# **•**TECAN•

Användarhandbok för



Dokumentnummer:	30225418
	2023-07
Dokumentutgåva:	1.5
Programvarurevisioner:	
magellan:	7.5
magellan Tracker:	7.5

CE



30225418 05



#### VARNING

#### LÄS NOGA IGENOM OCH FÖLJ INSTRUKTIONERNA I DENNA MANUAL INNAN DU ANVÄNDER INSTRUMENTET.

#### Anmärkning

Vi har gjort vårt bästa för att undvika fel i texter och diagram. Tecan Austria GmbH tar dock inte något ansvar för eventuella fel som kan förekomma i detta dokument.

Tecan Austria GmbH har som policy att förbättra produkter när ny teknik och nya komponenter blir tillgängliga. Tecan Austria GmbH förbehåller sig därför rätten att när som helst ändra specifikationerna baserat på adekvata verifieringar, valideringar och godkännanden.

Vi tar gärna emot eventuella kommentarer om innehållet i denna handbok.



#### Tillverkare

Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Österrike T: +43 6246 89 330 www.tecan.com E-post: office.austria@tecan.com

#### Upphovsrättsinformation

Innehållet i denna manual tillhör Tecan Austria GmbH och får inte kopieras, reproduceras eller överföras till annan person eller andra personer utan föregående skriftligt tillstånd.

Copyright © Tecan Austria GmbH Med ensamrätt. Tryckt i Österrike.

#### Deklaration om EC-godkännande

Tillhandahålls på begäran.

#### Avsedd användning för magellan

Se 1.2 Avsedd användning för magellan.

#### Om denna manual

magellan är ett universellt paket för reducering av data som används för att analysera data som genereras från analyser av mikroplattor. Detta är endast konstruerat för professionell användning.

Denna manual beskriver hur man:

- Installerar programvaran
- Använder programvaran

#### Anmärkning om skärmbilder

Det kan hända att versionsnumret som visas på skärmbilderna inte stämmer överens med numret på den aktuella utgåvan. Skärmbilder ersätts endast om innehållet som har att göra med programmet har ändrats.



#### Ändra historik för bruksanvisningar

Version	Ändringar	
V 1.2	Ny DoC (Declaration of Conformity, Konformitetsförklaring)	

#### Varningar, uppmaningar om försiktighet och anmärkningar

Följande typer av meddelanden används i den här publikationen. De belyser viktig information eller varnar användaren om eventuellt farliga situationer:





### Symboler

	Tillverkare
$\sim$	Tillverkningsdatum
CE	CE-märkning om överensstämmelse
ī	Se användarhandboken
IVD	Medicinteknisk produkt avsedd för in vitro-diagnostik
UK	Märkningen United Kingdom Conformity Assessed visar att produkten följer gällande bestämmelser i Storbritannien.
UDI	Unik enhetsidentifiering UDI-symbolen identifierar databäraren på etiketten.
REF	Katalognummer
SN	Serienummer
X	WEEE-symbol
	Kina RoHS-symbol



### Innehållsförteckning

1.	Inlee	dning		9
	1.1	Tilläm	npningsområde	9
	1.2	Avsed	dd användning för magellan	10
	1.3	Anvä	ndarprofil	10
		1.3.1	Professionell användare – administratörsnivå	10
		1.3.2	Slutanvändare eller rutinanvändare	10
	1.4	Speci	ifikationer	11
		1.4.1	Systemkrav	11
		1.4.2	Kompatibilitet med läsare	12
	1.5	Instal	lationsprocedur för programvara	12
		1.5.1	Automatisk installationsprogram	12
		1.5.2	Installation Qualification (Installationskvalificering) – IQ	13
		1.5.3	Operation Qualification (Driftskvalificering) – OQ	13
		1.5.4	Systemåterställning	14
		1.5.5	Automatisk avinstallation av programvara	15
		1.5.6	Cybersäkerhetsinformation	16
2.	Börj	ja arbe <sup>.</sup>	ta med magellan	17
	2.1	Anvä	ndargränssnitt – Guide-lista	17
	2.2	Komp	oonenter och termer – magellans grundläggande logik	19
		2.2.1	Filtyper som används med magellan	19
		2.2.2	Mapphantering	20
		2.2.3	Standardelement	22
		2.2.4	Knappen Help (Hjälp)	22
		2.2.5	Dialogrutan Välkommen	23
		2.2.6	Genvägslista	23
	2.3 Starta magellan		24	
		2.3.1	Start av standardversion	24
		2.3.2	Start av Trackerversion	24
	2.4	Anslu	ita ett instrument	25
		2.4.1	Ansluta till SUNRISE-instrument	25
		2.4.2	Ansluta till INFINITE F50-instrument	
		2.4.3	Licensiera magellan	
2	Inch	2.4.4	Registreringsguiden	28
3.	Inst	rument	tkontroll och installningar	
	3.1	Instru	imentalternativ	
		3.1.1	Movements (Förflyttning)	
		3.1.2	I emperature Control (I emperaturkontroll)	
	3.2	Instal	Iningar och servicealternativ	
		3.2.1	Change Instrument (Byt instrument)	
		3.2.2	Define Filter Slides (Definiera filternaliare)	
	3.3		les (Loggfiler)	
4.	Crea	ate/Edi	t a Method Wizard (Guiden Skapa/redigera en metod)	35
	4.1	Inledr	ning	35
	4.2	Defin	iera mätparametrarna	37
		4.2.1	Mättyper - SUNRISE	37
		4.2.2	Mätparametrar - SUNRISE	
		4.2.3	Mätparametrar - INFINITE F50	40
	4.3	Defin	iera utvärdering	42
		4.3.1	Fönstret Create/Edit Method Overview (Skapa/redigera metodöversikt)	
		4.3.2	Metodlayout: Hur man definierar en plattlayout	45



		4.3.3	Metodlayout: Koncentration, spädning och referensvärden	50
		4.3.4	Förhandsberäkning: Reducering av spektradata	53
		4.3.5	Transformed data (Transformerad data): Lägg till ny transformation	56
		4.3.6	Kinetik: Reducering av kinetiska data	59
		4.3.7	Kinetiska transformationer: Lägg till nya kinetiska transformationer	63
		4.3.8	Koncentrationer:Standardkurva	63
		4.3.9	Concentration transformation (Koncentrationstransformationer):	
			Lägg till nya koncentrationstransformationer	68
		4.3.10	Utvärdera data: Definition av cutoff	68
		4.3.11	Utvärdera data: QC-validering	70
		4.3.12	Datahantering: Dataexport	71
		4.3.13	Datahantering: Utskriven rapport	76
		4.3.14	Datahantering: Automatiserad datahantering	
		4.3.15	Miscellaneous (Ovrigt): Användarprompt	
		4.3.16	Miscellaneous (Ovrigt): Nummerformat	
		4.3.17	Miscellaneous (Ovrigt): Metodanteckningar	86
	4.4	Spara	metoden	86
		4.4.1	Lösenordsskydd av metoder	87
	4.5	Metod	er med flera plattor	
5.	Crea	te/Edit	a Sample ID List Wizard (Guiden Skapa/redigera en prov-ID-lista)	91
	5.1	Inledn	ing	91
	5.2	Create	e/Edit a Sample ID List (Skapa/redigera en prov-ID-lista)	
	0.2	5.2.1	Skapa ny prov-ID-lista)	
		5.2.2	Import/Edit a Sample ID List (Importera/redigera en prov-ID-lista)	
		5.2.3	Importera en prov-ID-lista	
		5.2.4	Spara prov-ID-listan	107
6.	Star	t Meası	urement Wizard (Guiden starta mätning)	109
•	6 1	Inledn		100
	6.0	Erbålla	n lý	140
	0.2			110
		6.2.1	Erhall radata med SUNRISE-Instrumentet	
	<u> </u>	0.2.2	Email radata med infiniti E F50-instrumentet	
	6.3	Use P	redefined Method (Anvand fordefinierad metod)	
	6.4	Start F	avorite (Starta Favorit)	111
	6.5	Starta	mätning med en fördefinierad eller favoritmetod	112
	6.6	Measu	irement Status (Mätstatus)	116
7.	Eval	uate Re	esults Wizard (Guiden Utvärdera resultat)	119
	7.1	Inledn	ing	119
	72	Select	a File (Väli en fil)	119
		721	Kriterier för filval	120
	73	Fönstr	et Workspace Overview (Översikt över arbetsvta)	121
	7.0	Elikon	Evoluete Deculte (Utvördere regultet)	100
	7.4		Monver	۲۲۵
		7.4.1		123
		7.4.2	Verktygsfältet: File (Fil)	124
		7.4.3	Verktygsfältet: Lott (Redigera)	120
		715	Fönstret nlattlavout	129 120
		74.0	Särskilda karaktärer	1∠9 120
		740	Kontrollfält Instrumentdata	130 120
		7 4 8	Kontrollfält Reducerade data	130
		1.7.0		100
		749	Kontrollfält Transformerad data	131
		7.4.9 7.4.10	Kontrollfält Transformerad data	131 131
		7.4.9 7.4.10 7.4.11	Kontrollfält Transformerad data Kontrollfält Kinetiska parametrar Kontrollfält Koncentrationer	131 131 132



		7.4.12 Kontrollfält Kvalitativa resultat	
		7.4.13 Kontrollfält Prov-ID:n	
		7.4.14 Kontrollfält Metodlayout	
		7.4.15 Kontrollfält QC-validering	136
		7.4.16 Kontrollfält Miscellaneous (Övrigt)	136
		7.4.17 Box för färgskala	137
		7.4.18 Sammanhangsberoende meny för en brunn	137
	7.5	Fliken Edit Method (Redigera metod)	142
	7.6	Spara de utvärderade resultaten	143
8.	Guid	len Bifoga underskrift	145
	8.1	Inledning	145
	8.2	Signera en fil	
9.	Ikon	en Miscellaneous (Övrigt)	
	91	Instrument control (Instrumentkontroll)	147
	9.1	File handling (Filhantering)	148
	5.2	9.2.1 Arkiv/filer	148
		9.2.2 Importera rådata	150
		9.2.3 Konvertera till	
		9.2.4 Konvertera från	
		9.2.5 Save Log Files (Spara loggfiler)	153
	9.3	Alternativ	
		9.3.1 Standarddatasökvägar	155
		9.3.2 Alternativ för kopiering/export	156
		9.3.3 Inställningar för plattvy	158
		9.3.4 Övrigt	159
	9.4	User Administration (Användaradministration) – magellan Tracker	161
		9.4.1 Lägg till/ändra användare (magellan Tracker)	
		9.4.2 Lägg till/ändra roll	
		9.4.3 Audit trail för användaradministration	
		9.4.4 Alternativ för användaradministration	
		9.4.5 Sammanfattning av användaradministration	
	9.5	User Administration (Användaradministration) - magellan Standard	169
		9.5.1 Lägg till/ändra användare (magellan Standard)	170
		9.5.2 Inloggning	170
		9.5.3 Andra användare	170
	9.6	Användarrättigheter	171
	9.7	Information om magellan	176
10.	Ytter	ligare funktioner för magellan Tracker	177
	10.1	User administration (Användaradministration)	177
		10.1.1 Audit Trail	177
	10.2	Filhantering	178
		10.2.1 Spara en fil	178
		10.2.2 Ändra en metodfil eller en arbetsyta	178
		10.2.3 Öppna en fil	178
		10.2.4 Öppna en fil som skapats på en annan PC – Lägg till HUID:er	178
	10.3	Systemets Audit Trail	179
11.	Berä	kningar	181
	11.1	Utvärdera resultaten – Beräkningsprocedur	
		11.1.1 Statistik	
	11.2	Reducering av spektradata	
		11.2.1 Matematisk beskrivning	



11.5 Exem 12.1 Ordlis ster märke	11.4.14 Beräkn 11.5.1 11.5.2 11.5.3 <b>pel på</b> Steg-fö 12.1.1 12.1.2 12.1.3 12.1.4 12.1.5 <b>sta me</b>	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> or-steg exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod) Kör metoden Evaluate the Result (Utvärdera resultat) Sammanfattning av definitionen av kvantitativ ELISA i magellan <b>d termer</b>	
11.5 Exem 12.1 Ordlis ster	11.4.14 Beräkn 11.5.1 11.5.2 11.5.3 <b>pel på</b> Steg-fö 12.1.1 12.1.2 12.1.3 12.1.4 12.1.5 <b>sta mee</b>	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> or-steg exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod) Kör metoden Evaluate the Result (Utvärdera resultat) Sammanfattning av definitionen av kvantitativ ELISA i magellan <b>d termer</b>	
11.5 Exem 12.1 Ordlis	11.4.14 Beräkn 11.5.1 11.5.2 11.5.3 <b>pel på</b> Steg-fö 12.1.1 12.1.2 12.1.3 12.1.4 12.1.5 <b>sta me</b>	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> prosets exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod) Kör metoden Evaluate the Result (Utvärdera resultat) Sammanfattning av definitionen av kvantitativ ELISA i magellan <b>d termer</b>	
11.5 <b>Exem</b> 12.1	11.4.14         Beräkn         11.5.1         11.5.2         11.5.3 <b>pel på</b> Steg-fö         12.1.1         12.1.2         12.1.3         12.1.4         12.1.5	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> or-steg exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod) Kör metoden Evaluate the Result (Utvärdera resultat) Sammanfattning av definitionen av kvantitativ ELISA i magellan	
11.5 <b>Exem</b> 12.1	11.4.14 Beräkn 11.5.1 11.5.2 11.5.3 <b>pel på</b> Steg-fö 12.1.1 12.1.2 12.1.3 12.1.4	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> pr-steg exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod) Kör metoden Evaluate the Result (Utvärdera resultat)	
11.5 <b>Exem</b> 12.1	11.4.14 Beräkn 11.5.1 11.5.2 11.5.3 <b>pel på</b> Steg-fö 12.1.1 12.1.2 12.1.3	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> or-steg exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod) Kör metoden	
11.5 <b>Exem</b> 12.1	Beräkn 11.5.1 11.5.2 11.5.3 <b>ipel på</b> Steg-fö 12.1.1 12.1.2	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> or-steg exempel: Kvantitativ ELISA Provsatsbeskrivning Create a Method (Skapa en metod)	
11.5 <b>Exem</b> 12.1	Beräkn 11.5.1 11.5.2 11.5.3 <b>pel på</b> Steg-fö 12.1.1	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b> or-steg exempel: Kvantitativ ELISA Provsatsbeskrivning	
11.5 <b>Exem</b> 12.1	Beräkn 11.5.1 11.5.2 11.5.3 <b>ipel på</b> Steg-fö	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) tillämpning or-steg exempel: Kvantitativ ELISA	
11.5 <b>Exem</b>	Beräkn 11.5.1 11.5.2 11.5.3 1 <b>pel på</b>	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden) <b>tillämpning</b>	205 205 205 206 206 <b>207</b>
11.5	11.4.14 Beräkn 11.5.1 11.5.2 11.5.3	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter) Calculation of IC Values (Beräkning av IC-värden)	205 205 205 206
11.5	11.4.14 Beräkn 11.5.1 11.5.2	ing i spädningsserien Upptäckt av spädningsserier Curve Parameter Calculation (Beräkning av kurvparameter)	205 
11.5	11.4.14 Beräkn 11.5.1	ing i spädningsserien Upptäckt av spädningsserier	205 205
11.5	Beräkn	ing i spädningsserien	205
	11.4.14		
		Viktning för fyra/fem parametrar – Marquardt/Polynomial Fit	204
	11.4.13	Fem parametrar – Marquardt	203
	11.4.12	Fyra parametrar – Marquardt	203
	11.4.11	Fyra parametrar	202
	11.4.10	LogitLog	200
	11.4.9	Akima	
	11.4.8	Kubisk spline	
	11.4.7	Polynomial (Polynom)	
	11.4.6	Non-Linear Regression (Icke-linjär regression)	
	11.4.5	l inear Regression (Liniär regression)	
	11 4 4	Point to point (Punkt-till-punkt)	190 196
	11 <u>4</u> 3	Felmeddelanden	190 106
	11.4.1	Parametrar för analystyn	
11.4		av anarys av Stanuarukurvur	
11 /	Typer c	Literinger	
	11.3.0		
	11.3.7	Andra runktioner	
	11.3.6	Eliminationstunktioner	
	11.3.5		
	11.3.4	Grundläggande funktioner	
	11.3.3	Formulelfunktioner	
	11.3.2	Formelvariabler	183
	11.3.1	Inledning	183
11.3	Hur ma	in skriver en formel	
	11.3	<ul> <li>11.3 Hur ma 11.3.1 11.3.2 11.3.3 11.3.4 11.3.5 11.3.6 11.3.7 11.3.8 11.3.9</li> <li>11.4 Typer a 11.4.1 11.4.2 11.4.3 11.4.4 11.4.5 11.4.6 11.4.7 11.4.8 11.4.9 11.4.10 11.4.11 11.4.12 11.4.13 11.4.14</li> </ul>	<ul> <li>11.3 Hur man skriver en formel.</li> <li>11.3.1 Inledning.</li> <li>11.3.2 Formelvariabler</li> <li>11.3.3 Formulelfunktioner</li> <li>11.3.4 Grundläggande funktioner.</li> <li>11.3.5 Statistiska funktioner</li> <li>11.3.6 Eliminationsfunktioner</li> <li>11.3.7 Andra funktioner</li> <li>11.3.8 Spektrafunctioner</li> <li>11.3.9 Exempel</li> <li>11.4 Typer av analys av standardkurvor</li> <li>11.4.1 Definitioner</li> <li>11.4.2 Parametrar för analystyp</li> <li>11.4.3 Felmeddelanden</li> <li>11.4.4 Point to point (Punkt-till-punkt)</li> <li>11.4.5 Linear Regression (Icke-linjär regression)</li> <li>11.4.6 Non-Linear Regression (Icke-linjär regression)</li> <li>11.4.7 Polynomial (Polynom)</li> <li>11.4.8 Kubisk spline</li> <li>11.4.9 Akima</li> <li>11.4.10 LogitLog</li> <li>11.4.11 Fyra parametrar – Marquardt</li> <li>11.4.14 Viktning för fyra/fem parametrar – Marquardt/Polynomial Fit</li> </ul>



### 1. Inledning

#### 1.1 Tillämpningsområde

magellan är en **programvara för styrning av läsare och dataanalys** för analysering av data som genereras av mätningar med en Tecan plattläsare i enkelläge för absorbans (t.ex. SUNRISE och INFINITE F50) för användning vid in vitro-diagnostik.



Anmärkning Det är viktigt att notera att enbart en korrekt installation av instrumentet och magellan-programvaran inte garanterar överensstämmelse med gällande förordningar. Motsvarande riktlinjer för processer och standardiserade tillvägagångssätt (SOP, standard operating procedures), inklusive validering och kvalitetskontroll måste också etableras.

**magellan** är utformad för att användas med de operativsystem och Microsoft Office-versioner som anges i kapitel 1.4.1 Systemkrav . Om ytterligare program installeras kan funktionaliteten i enlighet med IVD for Europa- och föreskrifterna i FDA inte garanteras.



Anmärkning För att förhindra missbruk av användarrättigheter och förfalskning av data rekommenderas användaradministratören att inte ha magellanrättigheter. Användaradministratören bör helst vara en del av ITavdelningen.



Kundernas systemadministratör är ansvarig för alla ändringar som görs i datorns operativsystem. Kunden måste se till att lämpliga användarrättigheter har ställts in för att undvika att data manipuleras eller raderas utanför magellan-programvaran.



Anmärkning

magellan är utformat för att användas med en dator. Det är inte avsett att integreras i ett nätverk.



Eventuella allvarliga incidenter som har inträffat i samband med enheten och programvaran ska rapporteras till tillverkaren och den behöriga myndigheten i den medlemsstat där användaren och/eller patienten är etablerad.

#### 1.2 Avsedd användning för magellan

**magellan** är en programvara för styrning av läsare och dataanalys för analysering av data som genereras av mätningar med en Tecan plattläsare i enkelläge för absorbans (t.ex. SUNRISE och INFINITE F50) för användning vid in vitro-diagnostik.

Programvaran **magellan** är avsedd att användas främst för absorbansmätningar och utvärdering av kvalitativa, semikvantitativa och kvantitativa IVD-analyser. Programvaran **magellan** kan programmeras av användaren; programvaran måste konfigureras av användaren i enlighet med analysinstruktionerna.

Produkten är avsedd för yrkesmässig laboratorieanvändning av utbildad personal. Produkten är inte avsedd för hemmabruk eller användning av privatpersoner.

magellan är tillgängligt i två versioner:

- magellan
- magellan Tracker

**magellan** Tracker erbjuder all nödvändig funktionalitet som krävs för att uppfylla föreskrifterna i FDA 21 CFR part 11 utöver **magellans** funktionalitet.



Anmärkning magellan är tillgängligt i två versioner: magellan och magellan Tracker.

För att underlätta hänvisningen i denna bruksanvisning är magellanversionen markerad som magellan Standard.

#### 1.3 Användarprofil

#### 1.3.1 Professionell användare – administratörsnivå

Administratören är en person med lämplig teknisk utbildning och motsvarade erfarenhet och färdigheter. Om produkten används som avsett kan denna person identifiera och undvika farliga situationer.

Administratören har omfattande kunskaper och kan instruera slutanvändare eller rutinanvändare gällande analysprotokoll med en Tecan-produkt inom dess avsedda användningsområde.

Vana vid datorprogram och goda kunskaper i engelska krävs.

#### 1.3.2 Slutanvändare eller rutinanvändare

Slutanvändaren eller rutinanvändaren är en person med lämplig teknisk utbildning och motsvarade färdigheter och erfarenhet. Om produkten används som avsett kan denna person identifiera och undvika farliga situationer.

Vana vid datorprogram och goda språkkunskaper i respektive nationellt språk vid installationsplatsen samt i engelska krävs.



Anmärkning Information om utbildningsdatum, utbildningarnas längd och hur ofta de genomförs finns tillgänglig via kundsupport. Adress och telefonnummer finns i användarhandboken och på nätet: http://www.tecan.com/customersupport

### 1.4 Specifikationer

#### 1.4.1 Systemkrav

Följande krav måste uppfyllas för att använda magellan:

	Stöds	Rekommenderat
PC	PC kompatibel med Windows med en processor kompatibel med Pentium på 1 GHz (Dual Core).	2 GHz (Dual Core)
Operativsystem	Windows 10 (32-bitars)	
	Windows 10 (64-bitars)	
	Versioner: Pro	
Minne	Windows 10 (32-bitars): 1 GB RAM	2 GB RAM
	Windows 10 (64-bitars): 2 GB RAM	4 GB RAM
Ledigt utrymme på hårddisk	3 GB	5 GB
Bildskärm	Super VGA-grafikkort	
Upplösning	1024 x 600 och högre	1920 x 1080
Färgdjup	256	
Mus	Microsoft-mus eller kompatibelt pekdon	
Kommunikation	1 x USB 2.0	2 x USB 2.0
		1 x RS232 (seriell)
Enheter	Windows 10:	
	DirectX 9 grafikkort med WDDM 1.0 eller högre	
.NET	Microsoft NET Framework 3.5: I Windows 10 uppmanas användaren att installera det nödvändiga .NET Framework (3.5) om det inte redan finns.	
Windows Installer	3.1	
	Om den här versionen inte finns kommer programmet för installation/uppgradering att installera den.	
Microsoft Excel	2007, 2010, 2013, 2016 (32-bitars),	2010 (32-bitars)
	2019 (32-bitars),	2019 (32-bitars)
	Microsoft Excel 365 (32-bitars)	
	Endast 32-bitarsutgåvor stöds!	
	Startutgåvor stöds INTE!	

#### 1.4.2 Kompatibilitet med läsare

Följande Tecan-läsare kan användas med magellan:

Instrumenttyper	Mätläge
SUNRISE	Absorbans
INFINITE F50	Absorbans

#### 1.5 Installationsprocedur för programvara

#### 1.5.1 Automatisk installationsprogram

magellan installeras av ett automatiskt installationsprogram. Alla nödvändiga komponenter installeras automatiskt.

Starta magellan\_Setup.exe för att påbörja installationen.

- I det första fönstret i guiden InstallShield Wizard visas de komponenter som krävs för att installera magellan. Klicka på Install (Installera) för att fortsätta.
   Windows 10: Om det nödvändiga .NET Framework (3.5) inte finns med, uppmanas användaren att starta installationen. Välj Install this feature (Installera den här funktionen) (en internetanslutning krävs). Om användaren hoppar över installationen kommer installationen av magellan inte att lyckas.
- Ett Välkomstfönster visas med information om programvaran och installationsprocessen. Klicka på Next (Nästa) för att fortsätta.
- Nästa fönster visar License Agreement (Licensavtalet). Läs igenom License Agreement (Licensavtalet) och klicka på I agree (Jag godkänner) för att acceptera och fortsätta.
- Därefter visas fönstret Customer Information (Kundinformation). Om du redan har köpt programvaran anger du ditt serienummer nu och klickar på Next (Nästa). Om du inte har ett serienummer än klickar du fortfarande på Next (Nästa) för att fortsätta installationen. Du kan registrera dig senare (se kapitel 2.4.3 Licensiera magellan).
- Därefter visas fönstret Destination Folder (Målmapp).
   Standardinställningssökvägen visas. Klicka på Browse (Bläddra) för att ändra standarddestinationen om så önskas. Klicka på Next (Nästa) för att fortsätta.
- Därefter visas fönstret Language Selection (Språkval). Välj önskat språk.
- I fönstret **Use For Regulated Environments (Används i reglerade miljöer)**, välj magellan **Standard** eller magellan **Tracker**, beroende på vilken version du har beställt, för att installera programvaran. Klicka på **Next (Nästa)** för att fortsätta.
- Installationsprogrammet är nu redo för installation. Klicka på **Next (Nästa)** för starta installationen.
- Klicka på **Finish (Slutför)** för att avsluta installationen och stänga installationsprogrammet.

Programvaran kan startas via Windows Start-menyn genom att välja magellan.

Installationsprogrammet upptäcker automatiskt en tidigare installation av **magellan**. Den gamla versionen måste avinstalleras innan du kan installera en ny magellan-version. Avinstallationen kan utföras automatiskt med hjälp av installationsguiden (se även kapitel1.5.5 Automatisk avinstallation av programvara).





i

Anmärkning Det är väldigt viktigt att personen som installerar programvaran har administratörsrättigheter på datorn.

Anmärkning Om en äldre version av magellan ska installeras måste den aktuella versionen avinstalleras helt innan den äldre versionen kan installeras från början.

#### 1.5.2 Installation Qualification (Installationskvalificering) – IQ

Det är rekommenderat att använda det automatiska installationskvalificeringsprogrammet för att kontrollera att installationen av **magellan** har lyckats.

Starta magellan IQ.exe från Windows start-menyn: **Start > Programs > Tecan >** magellan IQ eller direkt från filsystemet i den standardinställningsvägen som anges i installationsprogrammet (vanligtvis, **C:\Program Files\Tecan\magellan**).

Installationskvalificeringsprogrammet upptäcker automatiskt den installerade versionen av **magellan**. Klicka på **Check (Kontrollera)** för att börja installationskvalificeringen.

IQ-verktyget visar tillståndet för alla installerade komponenter. Det finns tre möjliga tillstånd: Successful (Lyckad), Warning (Varning) och Failed (=Error) (Misslyckad (=Fel)). Om eventuella fel rapporteras, kontakta din lokala kundtjänst för support.

Efter slutförd kontroll är det möjligt att skapa en rapport som innehåller den information som visas av IQ-verktyget. Förutom informationen i IQ-verktyget innehåller rapporten ett signaturfält så att rapporten kan sparas och skrivas ut i revisionsändamål. Klicka på knappen **Report (Rapport)** i: **File > Report (Fil > Rapport)** för att skapa en rapport. Rapporten kan sparas som en PDF-fil och andra filtyper.

Klicka på **Cancel (Avbryt)** eller **Exit (Avsluta)** för att stänga installationskvalificeringsprogrammet.



#### Anmärkning

Installationskvalificeringen bör upprepas varje gång magellan installeras, uppdateras till en nyare version, eller om det underliggande systemet uppdateras eller ändras.

#### 1.5.3 Operation Qualification (Driftskvalificering) – OQ

Utöver kontrollen av installationskvalificeringen rekommenderar Tecan också att **magellans** beräkningsförmåga testas. För det här syftet innehåller installationsmediet för **magellan** en arbetsyta och en rapportfil.

Öppna OQ-arbetsytan för den installerade **magellan**-versionen och skriv ut rapporten. Den nyligen utskrivna rapporten måste vara identisk med motsvarande OQ PDF-rapport (förutom tid och datum).

Om de två rapporterna inte är identiska, kontakta din lokala återförsäljare för support.



#### 1.5.4 Systemåterställning

För att reparera en skadad installation av **magellan** efter ett maskinvarufel är det viktigt att påbörja säkerhetskopieringenen av data i förväg.

#### Systemåterställning med säkerhetskopieringsprogram för data

Säkerhetskopiering av data utförs vanligtvis med säkerhetskopieringsprogram för data. Detta gör det möjligt att spara hela systemet, inklusive alla data, så att det kan återställas när som helst.

En detaljerad beskrivning och exakta instruktioner finns i Windows Help (för Windows säkerhetskopieringspaket) eller i den medföljande dokumentationen.

**magellan**-licensen är bunden till datorns maskinvara, vilket innebär att den kan bli ogiltig efter ett maskinvarubyte. Du bör därför kontrollera att din licens fortfarande är giltig efter en systemåterställning. Detta kan göras i dialogrutan **About magellan (Information om magellan)** genom att starta registreringsguiden och kontrollera om det registrerade serienumret och licensnumret fortfarande är giltiga. Om du får ett felmeddelande ska du kontakta Tecans kundsupport för att få ett nytt licensnummer.

#### Manuell systemåterställning

I de fall där en fullständig återställning av systemet inte är möjlig med säkerhetskopieringsprogrammet (t.ex. på grund av att maskinvaruförändringarna är för omfattande eller för att en annan version av Windows-operativsystemet installerades på den nya datorn), måste **magellan** installeras om och flera filer måste kopieras från en säkerhetskopia.

Då består systemåterställningsprocessen av följande steg:

- 1. Installera magellan
- 2. Starta magellan och definiera magellan-administratören
- 3. Ange användaren och användarrättigheterna på nytt.
- 4. Kopiera magellan.ini-filen från säkerhetskopian.
- 5. Kopiera magellan-metoderna från säkerhetskopian.
- 6. Kopiera prov-ID-listorna från säkerhetskopian.
- 7. Kopiera arbetsytorna från säkerhetskopian.
- 8. Kopiera exporterad data från säkerhetskopian.

När maskinvaruförändringarna och en omlicensering för magellan Tracker har slutförts får användaren information om att filerna kommer från en annan dator när användaren öppnar metodfilerna och arbetsytorna som skapats under den gamla licensen. För att lösa detta problem följer du den detaljerade beskrivningen i kapitel 10.2.4 Öppna en fil som skapats på en annan PC – Lägg till HUID:er.

#### Datasäkerhet

För att se till att stegen i systemåterställningen ovan kan slutföras snabbt bör en säkerhetskopia upprättas i tid. Säkerhetskopian bör innehålla **magellan**-metoder, arbetsytor, prov-ID-listor, exporterad data, audit trail-filer, system audit trail-filer, UserManagement.xml-filen, filerna i katalogen Pdfx och filen magellan.ini. Innehållet i säkerhetskopian ska alltid vara uppdaterat. Tecan rekommenderar att du använder alternativet **Auto Archiving (Automatisk arkivering)** för säkerhetskopiering av metoder, arbetsytor och prov-ID-listor. Du hittar information om Auto Archiving (Automatisk arkivering) i kapitel 9.2.1 Arkivfiler/Automatic Archiving (Automatisk arkivering).



magellan.ini	C:\Users\Public\Documents\Tecan\magellan
Metoder	Standardsökväg C:\Users\Public\Documents\Tecan\magellan\mth eller enligt inställningen i dialogrutan <b>Options (Alternativ)</b> .
Arbetsytor	C:\Users\Public\Documents\Tecan\magellan\wsp eller enligt inställningen i dialogrutan <b>Options (Alternativ)</b> .
Prov-ID-listor	C:\Users\Public\Documents\Tecan\magellan\smp eller enligt inställningen i dialogrutan <b>Options (Alternativ)</b> .
Exporterad data	C:\Users\Public\Documents\Tecan\magellan\asc eller enligt inställningen i dialogrutan <b>Options (Alternativ)</b> .
Audit trail-filer	Sökväg som visas i dialogrutan Audit Trail (under användaradministration).
Systemets Audit Trail	Sökväg enligt 10.3 Systemets Audit Trail
Användarhantering	C:\ProgramData\Tecan\Tecan User Management\v1.1\UserManagement.xml
Pdfx-filer	C:\Users\Public\Documents\Tecan\Pdfx

#### Sökvägar i magellan – Windows 10

#### 1.5.5 Automatisk avinstallation av programvara

Programvaran **magellan** kan tas bort med hjälp av Windows standardrutin för avinstallation:

- Stäng magellan
- Välj Add/Remove Programs (Lägg till/ta bort program) från Inställningar Kontrollpanelen i Windows start-meny.
- Välj ikonen för magellan och klicka på Remove (Ta bort).

Under borttagningen får du frågan om delade komponenter ska tas bort eller inte. Om **Yes (Ja)** väljs, kommer **magellan** att avinstalleras helt och andra Tecanprogram kommer inte längre att fungera. Det skadar inte ditt system om du låter dessa komponenter vara kvar. Om du är osäker på vad du ska göra föreslår vi att du inte tar bort de delade komponenterna. När du klickar på **Yes (Ja)** eller **No (Nej)** avinstalleras programvaran **magellan**.



#### Anmärkning

Vid borttagning av alla delade komponenter tas även data för användaradministration bort. Vissa Tecan-program, till exempel EVOware, kommer inte längre att fungera och måste installeras om.

#### 1.5.6 Cybersäkerhetsinformation

#### FÖRSIKTIGHET

#### DET ÄR ANVÄNDARENS ANSVAR ATT VIDTA NÖDVÄNDIGA FÖRSIKTIGHETSÅTGÄRDER FÖR ATT SKYDDA SYSTEMET FRÅN CYBERSÄKERHETSHOT.

VID INSTALLATION AV VIRUSSKYDDSPROGRAM ELLER SÄKERHETSRELATERADE UPPDATERINGAR AV OPERATIVSYSTEMET BÖR DU FÖLJA REKOMMENDATIONERNA FRÅN DEN LOKALA IT-AVDELNINGEN. SÄKERHETSKOPIERA MAGELLAN-FILER INNAN DU PÅBÖRJAR UPPDATERINGAR. KÖR INSTALLATIONSKVALIFICERING (IQ), DRIFTSKVALIFICERING (OQ) OCH PRESTANDAKVALIFICERING (PQ) AV MAGELLAN EFTER VARJE SYSTEMÄNDRING.

#### Säkerhetskopiera dina data

Vikten av att säkerhetskopiera filer som kan återhämtas i händelse av en ransomware-attack kan inte överdrivas. Se till att skapa destruktionsbeständiga säkerhetskopior av dina kritiska system och data. Det finns många verktyg och tjänster tillgängliga för säkerhetskopiering, återställning (av tidigare filversioner) och återställning av filer. Se också till att regelbundet testa att säkerhetskopiorna fungerar.

Använd magellans arkivfunktion eller egna verktyg för att säkerhetskopiera magellan-datafiler. För detaljerad beskrivning, se kapitel 9.2.1 Arkivfiler.

#### Hantera och kontrollera privilegierad tillgång till data

För att minimera risken för att autentiseringsuppgifter avslöjas och missbrukas bör du implementera unika lokala administratörslösenord på alla system, separera och skydda privilegierade konton och minska de omfattande behörigheterna i fildatabaserna.

Använd standad Windows-operatörer (inte administratörer) för att använda magellan.

Använd magellan Tracker-funktionen för att kontrollera åtkomst och ytterligare skydda programmet.

#### Driftskvalificering

Utöver kontrollen av installationskvalificeringen rekommenderar Tecan också att magellans beräkningsförmåga testas. För detaljerad beskrivning, se kapitel 1.5.3 Operation Qualification (Driftskvalificering) – OQ



#### Börja arbeta med magellan 2.

magellans standardguider representerar arbetsflödesmoduler. De är steg-försteg-guider för komplexa procedurer.

I vissa fall finns menyer tillgängliga i rubrikfältet. Menu (Meny) erbjuder ett konventionellt sätt att använda programvaran: det relevanta menyalternativet väljs från huvudmenyerna. Alla efterföljande åtgärder startas omedelbart, annars visas en dialogruta där ytterligare val eller poster kan göras.

#### 2.1 Användargränssnitt – Guide-lista

Den huvudsakliga typen av användargränssnittet i magellan är guiden.

När du har startat magellan visas Wizard List (Guide-listan):

$\cap$	What do you want to do?
Start measurement	<ul> <li>The Start Measurement wizard helps you to perform a measurement.</li> </ul>
Evaluate results	O You can either use a method or obtain raw data
Attach signature	0
Create/edit a sample ID list	•
Create/edit a method	•
/	

Varje guide kan startas antingen genom att dubbelklicka eller genom att välja den och klicka på knappen Next (Nästa).

#### Start Measurement Wizard (Guiden starta mätning)

Start Measurement wizard (Guiden Starta mätning) inkluderar följande alternativ:

- Obtain Raw Data (Erhåll rådata) används för att snabbt och enkelt generera rådata genom att ställa in de nödvändiga mätparametrarna och sätta igång en mätning.
- Use Predefined Method (Använd fördefinierad metod) används för att utföra mätningar baserat på metoder som skapats tidigare.
- Start favorite (Starta favorit) används för att välja en av de mest använda metoderna från listan med numrerade ikoner.

För detaljerad beskrivning, se kapitel 6 Start Measurement Wizard (Guiden starta mätning).

Efter slutförd mätning skapas en arbetsyta.

#### Evaluate Results Wizard (Guiden Utvärdera resultat)

**Evaluate Results wizard (Guiden Utvärdera resultat)** används för att visa rådata och för att utvärdera resultaten. Du kan se utvärderingsparametrarna och utvärdera data på nytt.

För detaljerad beskrivning, se kapitel 7 Evaluate Results Wizard (Guiden Utvärdera resultat).

#### Attach Signature Wizard (Guiden Bifoga underskrift)

Attach Signature wizard (Guiden Bifoga underskrift) används för att signera metoder och arbetsytor. Den här metoden är bara tillgänglig med magellan Tracker.

För detaljerad beskrivning, se kapitel 8 Guiden Bifoga underskrift.

#### Create/Edit a Sample ID List Wizard (Guiden Skapa/redigera en prov-ID-lista)

Create/Edit a Sample ID list wizard (Guiden Skapa/redigera en prov-ID-lista) används för att skapa nya och redigera befintliga prov-ID-listor.

För detaljerad beskrivning, se kapitel 5 Create/Edit a Sample ID List Wizard (Guiden Skapa/redigera en prov-ID-lista).

#### Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)

Create/edit a method wizard (Guiden Skapa/redigera en metod) används för att definiera eller redigera metoder.

För detaljerad beskrivning, se kapitel 4 Create/Edit a Method Wizard (Guiden Skapa/redigera en metod) .

#### Ikoner



#### Ikon: Ändra nuvarande användare

Om användaradministrationen är aktiv (se kapitel 9.4 User Administration (Användaradministration) – magellan Tracker och kapitel9.5 User Administration (Användaradministration) - magellan Standard) klickar du på den här ikonen för att logga ut den nuvarande användaren och logga in en ny användare.

Med **magellan** Standard är användaradministrationen valfri. Med **magellan** Standard är användaradministrationen obligatoriskt.



#### Ikon: Miscellaneous (Övrigt)

Klicka på ikonen Miscellaneous (Övrigt) för att välja följande alternativ:



Instrument control (Instrumentkontroll)



File handling (Filhantering)

g (Filhantening)

About **magellan** (Information om magellan)

Options (Alternativ)

User administration (Användaradministration)



- Instrument control (Instrumentkontroll) ger snabb tillgång till flera instrumentfunktioner, service och inställningsalternativ. Se avsnitt 3 Instrumentkontroll och inställningar..
- File handling (Filhantering) används för att flytta filer till en arkivplats, konvertera filer från eller till en annan magellan-version och för att importera rådata från en ASCII-fil. Se avsnitt 9.2 File handling (Filhantering).
- Options (Alternativ) används för att anpassa vissa standardinställningar som sökväg för skapade filer, alternativ för urklipp och Excel-kopiering, plattvyn och guiden övrigt, uppstart, inställningar för språk och lösenord. Se avsnitt 9.3 Alternativ.
- User administration (Användaradministration) används för att lägga till eller inaktivera användare och för att ställa in eller ändra användarrättigheter. Se kapitel 9.4 User Administration (Användaradministration) – magellan Tracker, respektive kapitel 9.5 User Administration (Användaradministration) - magellan Standard.
- About magellan (Information om magellan) tillhandahåller licensinformation och information om den installerade versionen och de installerade komponenterna. Registrering kan begäras genom att starta **Register wizard** (Registreringsguiden) (se kapitel 2.4.3 Licensiera magellan).

Stäng fönstret Miscellaneous (Övrigt) för att gå tillbaka till wizard list (guidelistan).

### Ikon: Temperature control (Temperaturkontroll)

Använd den här ikonen för att ställa in måltemperaturen för det anslutna instrumentet.

För detaljerad beskrivning, se kapitel 3.1.2 Temperature Control... (Temperaturkontroll)



#### Ikon: Move plate (Flytta platta)

Använd den här ikonen för att placera platthållaren i eller flytta den ut ur instrumentet.

#### 2.2 Komponenter och termer – magellans grundläggande logik

#### 2.2.1 Filtyper som används med magellan

Följande filtyper används med magellan.

**Som standard** lagras alla filtyper som är associerade med **magellan** i underkataloger i lämplig katalog:

...\All Users\Documents\Tecan\

#### 2. Börja arbeta med magellan

Underkatalogerna visas i tabellen nedan:

Type of File (Typ av fil)	File Extension (Filändelse)	Directory (Katalog)
Workspace (Arbetsyta)	.wsp	\magellan\wsp
Method (Metod)	.mth	\magellan\mth
Sample ID List (Prov-ID- lista)	.smp	\magellan\smp
Export Files (Exportfiler)	.asc	\magellan\asc
Standard Curve (Standardkurva)	.std	\magellan\wsp
Plate Definition (Plattdefinition)	.pdf .pdfx	\Reader\Pdf \Pdfx

För en detaljerad beskrivning av metoder, prov-ID-listor, arbetsytor och standardkurvor, se kapitel 4 Create/Edit a Method Wizard (Guiden Skapa/redigera en metod), 5.2.3 Importera en prov-ID-lista och 7 Evaluate Results Wizard (Guiden Utvärdera resultat).





Anmärkning INFINITE F50-instrument använder filformatet .pdfx, medan SUNRISEinstrument använder .pdf.

Anmärkning De tillgängliga menyerna och verktygsfältet varierar beroende på vilken typ av fil som för närvarande är öppen.

#### 2.2.2 Mapphantering

I **magellan**, är det möjligt att skapa en användarspecifik mappstruktur för att spara filer (metoder, arbetsytor, standardkurvor och prov-ID-listor) i vilken mapp som helst i Windows Explorer. Nya mappar kan skapas genom att högerklicka och välja **New folder (Ny mapp)** från den sammanhangsberoende menyn. Filer och mappar kan flyttas/klippas/kopieras på samma sätt som i Windows Explorer.

#### Symboler för filtyper

**magellan** använder följande symboler för metoder, arbetsytor, standardkurvor och prov-ID-listor i fönstret för filvyn:

2	Method (Metod)
	Standard Curve (Standardkurva)
(I)	Workspace (Arbetsyta)
	Password protected method (Lösenordsskyddad metod)
P	Sample ID List (Prov-ID-lista)



#### magellan Tracker

Standardsökvägar för att spara nyligen skapade filer kan ställas in via: Hemsidan för Wizard list (Guide-lista)  $\rightarrow$  knappen Miscellaneous (Övrigt)  $\rightarrow$  knappen Options (Alternativ) $\rightarrow$  fliken Paths (Sökvägar).

Användare kan skapa nya mappar i vilken mapp som helst i Windows Explorer under lagringsprocessen.

I fönstret **Save (Spara)** öppnas mappen som anges av filens standardsökväg automatiskt när en ny fil skapas. Om en befintlig fil ändras öppnas sökvägen till den aktuella platsen för filen. Det är dock möjligt att spara filer i vilken mapp som helst i Windows Explorer, eller i en nyligen skapad mapp.

#### magellan Tracker

Användare av **magellan** med administratörsrättigheter kan ställa in standardsökvägar för att spara nyligen skapade filer via: Hemsidan för **Wizard list (Guide-lista)**  $\rightarrow$  knappen **Miscellaneous (Övrigt)**  $\rightarrow$  knappen **Options** (Alternativ) $\rightarrow$  fliken **Paths (Sökvägar)**. Dessa standardsökvägar gäller för alla användare.

Användare kan skapa nya undermappar endast med Windowsadministratörsrättigheter och endast i den mapp som anges i standardsökvägen under lagringsprocessen.

I fönstret **Save (Spara)** öppnas mappen som anges av filens standardsökväg automatiskt när en ny fil skapas. Om en befintlig fil ändras öppnas sökvägen till den aktuella platsen för filen. Det är dock endast möjligt att spara filer i standardmappen, eller i en befintlig eller nyligen skapad mapp.



#### 2.2.3 Standardelement

Varje guide visar i tur och ordning ett antal fönster där all nödvändig information, inställningar och möjligheter till datainmatning finns.

#### Standardelement i en magellanguide

För att navigera från ett fönster till ett annat finns det flera knappar längst ner på guiden:

Knappen <b>Back</b> (Bakåt)	Knappen <b>Back</b> (Bakåt) används för att navigera tillbaka till det föregående fönstret inom en guide.
Knappen <b>Next</b> (Nästa)	Knappen <b>Next</b> (Nästa) används för att navigera framåt till nästa fönster inom en guide.
Knappen <b>Finish</b> Slutför)	Knappen <b>Finish</b> (Slutför) används för att navigera vidare till fönstret Save (Spara) i guiden.
Knappen <b>Save</b> (Spara)	Knappen <b>Save</b> (Spara) finns endast i det sista fönstret i en guide och ersätter knappen <b>Next</b> (Nästa). Den används för att avsluta guiden och spara alla ändringar, eller för att starta en process.
Knappen <b>Cancel</b> Avbryt)	Knappen <b>Cancel</b> (Avbryt) används för att avsluta en guide utan att spara några ändringar i inställningar eller dokument.
Knappen <b>Help</b> (Hjälp)	Knappen <b>Help</b> (Hjälp) öppnar hjälpfönstret.

#### Standardelement i Microsoft Windows

Knappen	Den här knappen bekräftar inställningarna, tillämpar och sparar
<b>OK</b>	ändringarna och stänger dialogrutan.
Knappen	Den här knappen stänger dialogrutan utan att spara några
<b>Cancel</b> Avbryt)	ändringar i inställningar eller dokument.
Knappen	Klicka på knappen <b>Help</b> (Hjälp) för att öppna <b>magellan</b>
<b>Help</b> (Hjälp)	onlinehjälp.

#### Information om statusfältet

Statusfältet visar följande information:

- Aktuell information om kommando
- Användarnamn på den inloggade användaren.
- Namn på det anslutna instrumentet. Till exempel: Sunrise
- Metod: mätläge och enhet. Till exempel: Absorbans [OD]
- Arbetsyta: datum och tid för mätningen. Till exempel: 27.11.2002 14:13:03 Antal utvalda brunnar. Till exempel: 3 brunnar utvalda
- Information om tangentbordets status: aktivitet f
  ör tangentbordets v
  äxelknappar: CAP (caps lock), NUM (Numeric block lock), SCRL (Scroll lock)
- Ikon för status för instrumentanslutning

#### 2.2.4 Knappen Help (Hjälp)

Klicka på knappen **Help (Hjälp)** eller tryck på "F1" för att öppna **magellan** onlinehjälp.



#### 2.2.5 Dialogrutan Välkommen

Varje guide börjar med dialogrutan **Welcome (Välkommen)**, som ger en kort introduktionsbeskrivning av den procedur som ska utföras.

Rensa kryssrutan **Show welcome page (Visa välkomstsida)** för att dölja välkomstsidor när du startar guider i framtiden.

#### 2.2.6 Genvägslista

SHIFT+B	Knappen Back (Bakåt)
SHIFT+N	Knappen Next (Nästa)
ESC	Knappen Cancel (Avbryt)
ENTER	Knappen Next (Nästa) eller Finish (Slutför), om det är ett aktivt fönster
F1	Hjälpmeny
CTRL+C eller CTRL+INSERT	Copy (Kopiera)
CTRL+V eller SHIFT+INSERT	Paste (Klistra in)
CTRL+X	Cut (Klipp ut)
CTRL+Y	Redo (Gör om)
CTRL+Z	Undo (Ångra)
DEL	Radera innehållet i den aktiva brunnen (redigera prov-ID, redigera formel)
CTRL+SHIFT	Visa formeln för den utvalda brunnen när transformationsresultatet visas (Guiden Evaluate Results (Utvärdera resultat)).

#### 2.3 Starta magellan

#### 2.3.1 Start av standardversion

Utför följande steg för att starta magellan:

- 1. Se till att ett instrument är anslutet eller välj ett simulerat instrument.
- 2. I Windows **Start**-menyn väljer du programgruppen **Tecan** och klickar på ikonen **magellan**.
- 3. magellan startar.



Anmärkning magellan kan köras med ett anslutet instrument eller i demoläge, där det simulerar ett instrument. Om du vill ansluta till ett instrument (se kapitel2.4 Ansluta ett instrument), sätt på instrumentet innan du startar magellan.



#### Anmärkning Innan du börjar arbeta med magellan rekommenderar vi att du läser versionsinformationen för magellan för att se om det finns några fel i programmet.

#### 2.3.2 Start av Trackerversion



Anmärkning När magellan Tracker används måste systemet för användaradministration anpassas av en utsedd administratör, som ansvarar för att skapa användarkonton och tilldelning av användarrättigheter. Administratören måste alltid utbildas av Tecan eller en organisation auktoriserad av Tecan.

#### Logga in för första gången med magellan Tracker

När **magellan Tracker** startas för första gången visas en dialogruta som informerar användaren om att en **User Administrator (Användaradministratör)** måste skapas först. Klicka på **OK** och dialogrutan **Create Administrator (Skapa administratör)** visas.

Fyll i textfälten och klicka på **OK** för att spara inställningarna. Minst en användaradministratör måste skapas.



FÖRSIKTIGHET OM ANVÄNDARADMINISTRATÖRENS LÖSENORD GLÖMS BORT KAN INSTÄLLNINGARNA FÖR ANVÄNDARADMINISTRATIONEN INTE ÄNDRAS OCH MAGELLAN MÅSTE INSTALLERAS OM HELT.

#### VI REKOMMENDERAR ATT MINST TVÅ ANVÄNDARADMINISTRATÖRER SKAPAS.

För alla ytterligare användardefinitioner i användaradministrationen (t.ex. användarrättigheter), se kapitel 9.4 User Administration (Användaradministration) – magellan Tracker, och kapitel 9.5 User Administration (Användaradministration) - magellan Standard.



#### Lösenord

#### Urspungliga lösenord (endast magellan Tracker)

När en användare loggar in för första gången måste lösenordet tilldelat av administratören ändras.

Efter det att användaren har angivit **User name (Användarnamn)** och **Password (Lösenord)** och klickat på **OK**, visas meddelandet "Your initial password is only valid once. You have to change the password!" (Ditt ursprungliga lösenord är bara giltigt en gång. Du måste byta lösenord!) Klicka på **OK**. Fönstret **Change Password (Byt lösenord)** tvingar användaren att ange ett nytt lösenord.

#### Utgånget lösenord (endast magellan Tracker)

När gilitghetstiden för lösenordet har gått ut uppmanas användaren att ange ett nytt lösenord.

Se kapitel 9.4.4 Alternativ för användaradministration.

#### Inloggning

När användaradministrationen har aktiverats visas dialogrutan Login (Inloggning) varje gång magellan startar.

Dialogrutan Login (Inloggning) innehåller följande element:

Textrutan <b>Username</b> (Användarnamn)	Ange ditt Användar-ID
Textrutan Password (Lösenord)	Ange ditt Lösenord

Klicka på knappen GO (Gå) för att starta eller på knappen Cancel (Avbryt) för att avsluta magellan.

#### Programmet låst

Om programmet inte har använts under den angivna maximala tidsperioden (användardefinierad endast i **magellan** Tracker), kommer det att låsas för användning. Ett lösenord måste anges för att låsa upp programmet.

#### 2.4 Ansluta ett instrument

#### 2.4.1 Ansluta till SUNRISE-instrument

Anslut instrumentet till datorn enligt beskrivningen i bruksanvisningen för instrumentet och slå på instrumentet innan du börjar **magellan**.

#### Första gången magellan startas

Första gången magellan startas visas dialogrutan Setup Port (Portinställning):

Instrument	W		OK
Select:	SUNRISE	~	-
◯ Find any	10 10		Cancel
Port			Help
◯ Select:	Find any	~	
Find any			
O Demo mode			



Instrument	Välj det anslutna instrumentet i rullgardinsmenyn eller välj <b>Find any</b> (Leta upp).
Port	Välj lämplig <b>COM port</b> från rullgardinsmenyn eller välj <b>Find any</b> (Leta upp) och klicka på <b>OK</b> .

#### **Change Instrument (Byt instrument)**

Klicka på ikonen **miscellaneous (Övrigt)** i guide-listan och **instrument control** (Instrumentkontroll) i listan miscellaneous (Övrigt). Välj Change instrument (Byt instrument) från Setup & Service (Inställning och service) för att ansluta ett instrument eller för att ändra det anslutna instrumentet.

#### Demoläge

Om **inget instrument** är anslutet, välj det instrument som ska simuleras från rullgardinsmenyn i grupprutan **Instrument** och välj **Demo mode (Demoläge)** i grupprutan **Port**. Klicka på **OK**. **magellan** är nu i demoläge.

etup Port			×
Instrument	Terrane and the second s		OK
Select:	SUNRISE	~	C 1
◯ Find any			Lancel
Port			Help
O Select:	Find any	~	
◯ Find any			
Demo mode			

**Demo mode allowed (Demoläge tillåtet)** (i dialogrutan **Options (Alternativ)**) är valt som standard.

I demoläget är det möjligt att utföra alla funktioner förutom faktiska mätningar.

#### 2.4.2 Ansluta till INFINITE F50-instrument

#### Första gången magellan startas

Första gången magellan startas visas dialogrutan Setup Port (Portinställning):

Instrument			OK
Select:	infinite F50	~	
◯ Find any			Cancel
Port			Help
◯ Select:	Find any	~	
Find any			

Välj INFINITE F50 i grupprutan Instrument i dialogrutan Setup Port (Portinställning). Klicka på OK för att visa dialogrutan Connect to Instrument (Anslut till instrument).



La constanti Norma	<b>T</b>	01:	D-4
	Туре	Alias	FOR
ntinite FSU	READER	1601003620	USBU
dditionally connect to:			
Instrument Name			Port
] Show simulated instrument	s		

Välj instrumentet och klicka på OK.

#### Demoläge

Om **inget instrument** är anslutet, välj det instrument som ska simuleras från rullgardinsmenyn i grupprutan **Instrument** i dialogrutan **Setup Port (Portinställning)**:

Linter and Marian	<b>T</b> 222	Alizz	D-4	-
infinite F50	Reader	Simulation	SUSIM:	
dditionally connect to:				
Instrument Name			Port	
ConnectSimulator			Connect	
				_
Show simulated instruments		F50_4_FILTERS		

I demoläget markerar du kryssrutan **Show simulated instruments (Visa simulerade instrument)** och det önskade instrumentet i rullgardinsmenyn, klicka på instrumentets namn och klicka på **OK**.



#### 2.4.3 Licensiera magellan



Anmärkning

magellan kan användas oregistrerad under 30 arbetsdagar. Om magellan inte har registrerats efter denna period, kommer alternativen spara och utskrift att avaktiveras.

Licensieringen av **magellan** görs med hjälp av guiden **Register magellan** (**Registrera magellan**). Om **magellan** inte har blivit registrerad, kommer denna guide att automatiskt starta varje gång programmet startas.

Guiden Register magellan (Registrera magellan) kan också öppnas via ikonen miscellaneous (Övrigt) på startsidan för guide-listan. (miscellaneous  $\rightarrow$  About magellan  $\rightarrow$  Register).

För att kunna registrera **magellan**, behövs ett serienummer från magellan, maskinvarans (PC) identifikationsnummer (HUID) och ett licensnummer.

För att licensiera magellan finns det två möjligheter:

1. Programvaran har köpts. Serienumret levereras tillsammans med programvaran.

När serienumret har angetts i dialogrutan register **magellan** (Registrera magellan) fyller du i registreringsformuläret för att påbörja begäran om licensnummer. Det ifyllda registreringsformuläret skickas tillsammans med det individuella HUIDnumret till Tecan Austria. HUID-numret skapas av programvaran och läggs automatiskt till i registreringsformuläret. Det är relaterat till Windows systemenhetsnummer.

När du har fått licensnumret måste du starta registreringsguiden igen och ange numret. Guiden **Register magellan (Registrera magellan)** bekräftar licensnumret och visar en summering av användarinformationen: Klicka på knappen **Finish (Slutför)** för att slutföra registreringen. **magellans** funktioner kommer därefter att vara fullt åtkomliga för användaren.

 Programvaran levererades som en demoversion tillsammans med instrumentet. En registrering är möjlig efteråt (inget serienummer levereras).

För att köpa **magellan** och få ett serienummer väljer du alternativet **Order magellan to obtain a valid license (Beställ magellan för att erhålla en giltig licens)** och fortsätt med registreringen enligt beskrivningen ovan.



#### Anmärkning

Det är väldigt viktigt att personen som registrerar programvaran också har administratörsrättigheter för datorns operativsystem

#### 2.4.4 Registreringsguiden

Registreringsguiden börjar med dialogrutan **Welcome (Välkommen)**, som innehåller en kort beskrivning av guiden. Klicka på **Next (Nästa)** och dialogrutan **Serial Number (Serienummer)** visas.



#### Serienummer och licensnummer

#### Beställ ett serienummer för magellan

För att köpa **magellan** och få ett serienummer väljer du alternativet **Order magellan to obtain a valid license (Beställ magellan för att erhålla en giltig licens)**. Fönstret **Registration Form (Registreringsformulär)** visas då (se kapitelRegistreringsformulär nedan).

Register Magellan	×
Enter your Magellan serial numbe	r: -
	< Back Next > Cancel

#### magellan har redan köpts

Välj Enter your magellan serial number (Ange ditt serienummer från magellan). Ange serienumret som finns på magellans förpackning.

Efter att du har angett serienumret, klicka på **Next (Nästa)** och dialogrutan **License Number (Licensnummer)** visas.

gister magenan				
Enter your Magellan licens	e number:			
O Request a Magellan licens	se number because y	ou did not rece	aive one.	

Användare som installerar programvaran för första gången har ännu inget licensnummer och måste välja **Request a magellan license number because** you did not receive one (Begär ett licensnummer för magellan eftersom du ännu inte har fått något). Klicka på Next (Nästa) och fönstret **Registration** Form (Registreringsformulär) visas.



#### Registreringsformulär

Fyll i registreringsformuläret . Fälten som är markerade med en asterisk är obligatoriska.

Full Name:	Test User				
Company / Institution:	Tecan Austria				
Street Address 1:	Untersbergstrasse 1a				•
Street Address 2:		10			
ZIP, City / Town:	5082 Groedig	* Country:	Austria	~	
Telephone:		Fax:			
E-Mail:	test.user@tecan.com	-			
Area of Work (Life Scienc	e/Clinical Diagnostic/Othe	·):		~	
Used Tecan Instruments:	SUNRISE			Add	
Detection Modes	Absorbance		Add		

Klicka på **Next (Nästa)** och fönstret med licensavtalet visas. Läs den noggrant och klicka på knappen **I agree (Jag godkänner)**.

Klicka på Next (Nästa) och registreringsinformationen visas.

Your Registration Re	equest Info		
USER ADDRESS	Test User / Tecan A Untersbergstrasse la	Austria a	^
	5082 Groedia / Aust	tria	
APPLICATION:	Magellan, 7, 2, 1,	5	
WORKSTATION	94086-30335FB		~
<			>
E-Mail Form	Print Form	View Form	

Klicka på **E-Mail Form... (Skicka formulär via e-post)** för att skicka informationen med hjälp av standardprogrammet för e-post på din dator eller klicka på **Print Form... (Skriv ut formulär)** för att skriva ut registreringsformuläret för att sedan faxa det eller skicka det med vanlig post till Tecan. Klicka på **View Form... (Visa formulär)** för att visa registreringsformuläret i WordPad eller Notepad. Inom 24 timmar får du licensnumret.



#### Slutför licensiering av magellan

När du har fått Tecans licensnummer upprepar du stegen ovan och lägger till licensnumret. Klicka på **Next (Nästa)**. Registreringsguiden bekräftar licensnumret och visar en summering av användarinformationen.

Klicka på knappen **Finish (Slutför)**. Guiden **Register magellan (Registrera magellan)** stängs och **magellans** funktioner kommer att vara fullt åtkomliga för användaren.

### 3. Instrumentkontroll och inställningar.

Klicka på knappen Instrument Control (Instrumentkontroll) i dialogrutan Wizard List (Guide-lista) och dialogrutan Instrument Control (Instrumentkontroll) visas.

Beroende på vilket instrument som är anslutet aktiveras eller inaktiveras olika instrument-, inställnings- och servicealternativ.

#### 3.1 Instrumentalternativ

#### 3.1.1 Movements... (Förflyttning)

Detta öppnar dialogrutan **Movements (Förflyttning)** där det är möjligt att styra förflyttningar på platthållaren och filterhållaren.

#### 3.1.2 Temperature Control... (Temperaturkontroll)

Endast tillgänglig för instrument som är utrustade med temperaturkontroll.

Med detta alternativ kan användaren fastställa temperaturen i läsaren.

Den här dialogrutan är också tillgänglig via dialogrutan Start Measurement (Starta mätning) innan en mätning startas. (Se kapitel6.5 Starta mätning med en fördefinierad eller favoritmetod).

Temperature Control	
Current temperature Current temperature: 34.9 °C Refresh	OK Cancel Help
Target temperature  Temperature control on  Target temperature: 37 °C  Apply	Plate in

<b>Current</b>	Den aktuella temperaturen visas i motsvarande fält. Klicka
temperature	på knappen <b>Refresh</b> (Uppdatera) för att uppdatera den
Aktuell temperatur)	aktuella temperaturen.
<b>Target</b>	För att starta eller stoppa temperaturkontrollen markerar
temperature	eller avmarkerar du kryssrutan och anger måltemperaturen.
Måltemperatur)	Klicka på <b>Apply</b> (Tillämpa)
	för att skicka den angivna temperaturen till instrumentet och klicka slutligen på <b>OK</b> för att stänga dialogrutan.

#### 3.2 Inställningar och servicealternativ

#### 3.2.1 Change Instrument... (Byt instrument)

l rutan Setup & Service (Inställningar och service) klickar du på Change instrument... (Byt instrument) för att öppna dialogrutan Setup Port (Portinställning).

Detta alternativ gör det möjligt att ansluta **magellan** till ett instrument. Se kapitel2.4 Ansluta ett instrument för mer detaljer.



Anmärkning Om du ansluter ett annat instrument till datorn eller om du ändrar gränssnittsparametrarna måste du alltid välja det här menyalternativet.

Inställningarna används automatiskt nästa gång programmet startas.

#### 3.2.2 Define Filter Slides... (Definiera filterhållare)

Det går att definiera standardfilterhållare och anpassade filterhållare. Se motsvarande kapitel i instrumentets användarhandbok.

#### 3.3 Log Files (Loggfiler)

När du arbetar med magellan skapas loggfiler. Kommunikationen mellan programvaran och instrumentet och kommunikationen mellan programvarans komponenter lagras i dessa loggfiler.

De finns i följande sökvägar:

Loggfiler för magellans (kommunikation mellan enstaka komponenter i magellan

Windows 10:

C:\Users\Public\Documents\Tecan\LogFiles\

- Insamling av nödvändiga loggfiler
   Loggfiler kan sparas som zip-arkiv genom att välja knappen Save Logfiles...
   (Spara loggfiler) i dialogrutan File handling (Filhantering) (Miscellaneous
   → File Handling). Zip-arkivet kan nu namnges och sparas i en angiven katalog. I händelse av mät- eller statusfel i magellan, innehåller detta arkiv alla data om brunnar, status (t.ex. överflöde, "lamp low") eller beräkningsfel och kan enkelt skickas till din lokala kundtjänst för support. För mer information se kapitel 9.2 File handling (Filhantering) Spara loggfiler.
- Loggfiler för INFINITE F50 (kommunikation mellan magellan och INFINITE F50-instrumentet): Windows 10: C:\Lisers\Public\Documents\Tecan\LogEiles\magellan\V x v\Instrument Serial

 Loggfiler för SUNRISE (RdrOle.log; kommunikation mellan magellan och SUNRISE-instrumentet): Windows 10: C:\Users\CurrentUser\AppData\Local\Temp



## 4. Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)

#### 4.1 Inledning

Create/Edit a Method Wizard (Guiden Skapa/redigera en metod) används för att

- Skapa eller redigera metoder,
- Ställa in parametrar för mätning och utvärdering,
- Definiera plattlayouten,
- Välja format för den utskrivna rapporten och
- Ställa in parametrarna för automatiserad datahantering.

#### Sammanfattning av arbetsflöde

Klicka på Continue (Fortsätt) på Välkomstsidan av Create/edit a method wizard (Guiden Skapa/redigera en metod) I nästa fönster, välj

- Create new (Skapa ny) om du vill definiera en ny metod eller
- Edit (Redigera) om du vill ändra en befintlig metod.

I fönstret **Define Measurement Parameters (Definiera mätparametrar)** kan mätparametrar ställas in. Klicka på **Define Evaluation (Definiera utvärdering)** för att definiera plattlayouten, utvärderingsparametrar, utskrift och parametrar för automatiserad datahantering. I slutet av guiden sparas den nya eller ändrade metoden som en .mth-fil.



Anmärkning För en steg-för-steg-beskrivning om hur man skapar en metod, se Kapitel 12 Exempel på tillämpning.

Sida för filval

I guide-listan, klicka på knappen Create/Edit a method (Skapa/redigera en metod). Klicka på Next (Nästa) på Välkomstsidan av Create/edit a method wizard (Guiden Skapa/redigera en metod) och sidan File Selection (Filval) med följande element visas:

Knappen <b>Create new</b> (Skapa ny)	Välj knappen <b>Create New</b> (Skapa ny) för att skapa en ny metod.
Knappen <b>Edit</b> (Redigera)	Knappen <b>Edit</b> (Redigera) måste väljas för att redigera en befintlig metod.



Show (Visa) kombinationsruta	<ul> <li>I kombinationsrutan Show (Visa) kan du definiera ett filfilter för att endast få fram de metoder i listan som uppfyller de valda kriterierna:</li> <li>All files (Alla filer)</li> <li>Files from this instrument (Filer från det här instrumentet)</li> <li>My files (Mina filer): Det här alternativet är tillgängligt om användaradministrationen är aktiverad (den är alltid aktiverad i magellan Tracker).</li> <li>Signed files (Signerade filer): endast tillgängliga för magellan Tracker</li> <li>Example files (Exempelfiler): endast tillgängliga om de har installerats.</li> </ul>
Listan <b>ilename</b> (Filnamn)	Välj den metod som ska redigeras från listan <b>Filename</b> (Filnamn). Fältet <b>Remarks</b> (Kommentarer) bredvid varje filnamn innehåller en kort beskrivning av metoden, om den anges. Alla metoder som finns tillgängliga i standardmetodkatalogen visas (se kapitel9.3 Alternativ).
Knappen <b>Print Preview</b> (Förhandsgranskning av utskrift)	Klicka på knappen <b>Print Preview</b> (Förhandsgranskning av utskrift) för att öppna <b>dialogrutan Print Preview</b> (Förhandsgranskning av utskrift) där en förhandsgranskning av inställningarna för den valda metoden visas och en utskrift kan startas.



#### Anmärkning

Om ett SUNRISE-instrument är anslutet, men en INFINITE F50-metod är vald, visas dialogrutan Instrument Mismatch (Missmatchade instrument):

• Anslut till rätt instrument.


# 4.2 Definiera mätparametrarna

I fönstret **Measurement Parameters (Mätparameterar)** är det möjligt att ställa in alla nödvändiga parametrar för mätningen, inklusive mätläge, våglängder, läsningsläge, temperatur osv., beroende på typen av instrument som är ansluten.

När du är ansluten till ett INFINITE F50-instrument, se användarhandboken för icontrol för mer information om hur du definierar mätparametrar.

# 4.2.1 Mättyper - SUNRISE

Klicka på Make your selection (Gör ditt val) och dialogrutan Measurement type (Mättyp) visas:

Create/Edit a Method	×
	Measurement type
	Endpoint measurement     Well kinetic measurement     Well kinetic measurement
	Measurement parameters
Help Cancel <<< Back	DEFINE EVALUATION

Mättypen ska väljas från en lista med alternativknappar.

Fönstret **Measurement type (Mättyp)** innehåller följande element (beroende på instrumentet som är anslutet):

Grupprutan <b>Measurement type</b> (Mättyp)	<ul> <li>Välj Endpoint measurement (Slutpunktsmätningar) för att utföra en enstaka mätning.</li> </ul>
	<ul> <li>Välj Kinetic measurement (Kinetisk mätning) för att utföra kinetiska mätningar per platta med en angiven intervalltid.</li> </ul>
	<ul> <li>Välj Multilabel measurement (Mätningar med multipla inmärkningar) för att utföra mätningar med multipla inmärkningar, med olika mätparametrar.</li> </ul>

Klicka på Measurement parameters... (Mätparametrar) för att öppna dialogrutan Measurement Parameter (Mätparameter).

# 4.2.2 Mätparametrar - SUNRISE

Välj önskad **Measurement type (Mättyp)** och klicka på **Measurement parameters... (Mätparameterar)** för att öppna dialogrutan Measurement Parameter (Mätningsparameter) för justering av

- Barcode (Streckkod),
- Wavelengths (Våglängder),
- Measurement parameters (read mode) (Mätparameterar läsningsläge),
- Kinetics (Kinetik)
- Temperature (Temperatur) och
- Shaking (Skakning).

Klicka på knappen **OK** för att spara ändringarna och på knappen **Cancel (Avbryt)** för att avvisa dem.

Beroende på vilket instrument som är anslutet och mättypen som valts varierar dialogrutan **Measurement Parameters (Mätparameterar)** och de respektive flikarna som är tillgängliga.

Vid anslutning till ett Infinite-instrument, se användarhandboken för i-control.

#### Multilabel Measurement (Mätningar med multipla inmärkningar)

Om mättypen Multilabel measurement (Mätningar med multipla inmärkningar) har valts visas dialogrutan Multilabel Measurement (Mätningar med multipla inmärkningar)

För att skapa en lista med parameterfönster måste du klicka på knappen **New** (Ny). Dialogrutan **Measurement Parameters (Mätparameterar)** visas där ett nytt set mätparametrar kan definieras. Minst två set av mätparametrar måste skapas i dialogrutan multipla inmärkningar.



Anmärkning

Om Move plate out after measurement (Flytta ut plattan efter mätning) väljs när metoden körs, visar magellan en meddelanderuta där vätskorna i plattan kan ändras och mätningen kan fortsätta. I annat fall utförs mätningen utan paus.

Mätparametern **Comments (Kommentar)** kan användas för att märka mängden mätparametrar i listan.



Multilabel Measurement X New... No. Label Measurement parameters 
 I
 DEF-1
 Measurement mode: Absorbance;Measurem

 I
 2
 DEF-2
 Measurement mode: Absorbance;Measurem

 I
 3
 DEF-3
 Measurement mode: Absorbance;Measurem
 Properties. Delete < > Kinetic 2 Number of cycles: 000:00:00 hh:mm:ss Interval: Use minimum interval Cancel Help OK

Listan <b>Label (Märkning)</b>	I listan <b>Multilabel (Multipla inmärkningar)</b> listas de befintliga definitionerna av mätparametrar rad för rad. I kolumnen <b>No. (Nr.)</b> räknar ett referensnummer upp de befintliga definitionerna av mätparametrar och en liten ikon visas. Kolumnen <b>Label (Märkning)</b> visar ett anpassningsbart namn (i <b>Measurement Parameters –</b> <b>General Tab – Comment to this measurement</b> ). Kolumnen <b>Measurement parameters (Mätparameterar)</b> visar en sammanfattning av de valda mätparametrarna.	
Knappen <b>New (Ny)</b>	Klicka på knappen <b>New (Ny)</b> och mätparametrarna för varje ny mätning kan definieras. Dialogrutan <b>Measurement Parameters (Mätparametrar)</b> visas. Definiera namn för varje märkning. I dialogrutan <b>Measurement Parameters (Mätparameterar)</b> används namnet som skrivits i kommentarsfältet på fliken <b>General</b> (Allmänt) som märkning. Alla parametrar måste anges för att definiera en ny mätning.	
Knappen <b>Properties</b> (Egenskaper)	Klicka på knappen <b>Properties (Egenskaper)</b> och alla <b>measurement parameters (Mätparameterar)</b> för den markerade märkningen i listan över mätparametrar visas och kan redigeras.	
Knappen <b>Delete (Radera)</b>	Klicka på <b>Delete (Radera)</b> eller tryck på <b>DEL</b> för att radera den/de markerade mätparametern/mätparametrarna från fältet med multipla inmärkningar.	
Number of cycles (Antal cykler)	Definierar antalet kinetikcykler för en kinetisk mätning med multipla inmärkningar.	
Interval (Intervall)	Definierar det kinetiska intervallet eller pausen mellan märkningarna (om endast en cykel definieras).	

Dialogrutan Multilabel (Multipla inmärkningar) innehåller följande element:

# 4.2.3 Mätparametrar - INFINITE F50

#### Definiera slutpunktsmätningar

Följande exempel beskriver en **Absorbance Endpoint Measurement** (Slutpunktsmätning för absorbans) i alla brunnar i en mikroplatta med 96 brunnar.

- 1. Välj en mikroplatta med 96 brunnar från rullgardinsmenyn **Plate definition** (Plattdefinition).
- 2. Som standard väljs alla brunnar i mikroplattan med 96 brunnar för mätning.
- 3. Ange önskad mätnings- och referensvåglängd.

		2
	tals.	
ŏŏŏŏŏŏŏŏŏŏŏ		
Absorbance		3
V Absorbance	Label	3
	Label Name: Labelt v	3

## Definiera mätningar med multipla inmärkningar

Följande exempel beskriver en **Absorbance Multilabel Measurement (Mätning med multipla inmärkningar för absorbans)** i alla brunnar i en mikroplatta med 96 brunnar. Tre absorbansmärkningar ska mätas.

- 1. Välj en mikroplatta med 96 brunnar från rullgardinsmenyn **Plate definition** (Plattdefinition).
- 2. Som standard väljs alla brunnar i mikroplattan med 96 brunnar för mätning.
- 3. Ange önskad mätningsvåglängd.
- 4. Infoga 2 eller fler **Absorbance (Absorbans)**-element och ange mätningsvåglängderna.

✓ Part of Plate		
	Datak	
Absorbance		
Wavelength	Label	
Measurement: 405 nm ~	Name: Label1 ~	
Reference: 405 rm		
Absorbance		
Wavelength	Label	
Measurement: 492 nm 🗸 🗸	Name: Label2 ~	
Reference: 405 nm		
🖞 🔻 Absorbance		
Absorbance Wavelength	Label	
Vavelength Measurement 620 nm ~	Label Name: Label3 ✓	



#### Definiera kinetiska mätningar

Följande exempel beskriver en kinetisk mätning i en mikroplatta med 96 brunnar.

- 1. Välj en mikroplatta med 96 brunnar från rullgardinsmenyn **Plate definition** (Plattdefinition).
- 2. Infoga ett **Kinetic Cycle (Kinetikcykel)**-programelement mellan delen av plattan och absorbans-elementet.
- 3. Cycles (Cykler)/Number of cycles (Antal cykler): 50
- 4. Kinetic interval (interval between measurements) (Kinetikintervall (intervall mellan mätningar)): välj **Use kinetic interval (Använd kinetikintervall)** och ange: 2 minuter 30 sekunder
- 5. Definiera **Absorbance (Absorbans)**-elementet genom att ange önskad mätningsvåglängd.

🐐 🔻 Part of Plate		2
$\begin{array}{c ccccccccccccccccccccccccccccccccccc$	Details	
😡 🔻 Kinetic Cycle		3
Cycles	Kinetic Interval	
Number of cycles: 50 +	Use kinetic interval:	
O Duration	Time: 00:02:30   ⊕  (hh:mm:ss)	
	O Time: 150000.2 ma	
Absorbance		4
Wavelength	Label	
Measurement: 492 nm	V Name: Label1 V	
Reference: 405 nm		

# 4.3 Definiera utvärdering

# 4.3.1 Fönstret Create/Edit Method Overview (Skapa/redigera metodöversikt)

I fönstret **Define Evaluation (Definiera utvärdering)** definierar användaren plattlayouten, transformationer och beräkningar, väljer format för den utskrivna rapporten och ställer in parametrarna för automatiserad datahantering.



# Verktygsfältet

Överst i fönstret visas ett **Verktygsfält** med de vanligaste funktionerna beroende på de aktuellt valda alternativen:

Knappen <b>Undo</b> (Ångra)	Klicka på <b>Undo</b> (Ångra) för att ångra en tidigare åtgärd.
Knappen <b>Redo</b> (Gör om)	Klicka på <b>Redo</b> (Gör om) för att göra om den ångrade åtgärden.
Knappen <b>Select all unused</b> (Välj alla oanvända)	Endast i läget plattlayout: Alla oanvända brunnar på plattan är markerade.
Knappen <b>Zoom to 10%</b> (Zooma till 10 %)	Detta ställer in plattlayoutvyn på 10 %.
Knappen <b>Zoom to 100%</b> (Zooma till 100 %)	Detta ställer in plattlayoutvyn på 100 %.
Knappen <b>Zoom mode</b> (Zoomläge)	Använd knappen <b>Zoom mode</b> (Zoomläge) för att zooma in i det markerade området. Om den är aktiverad kan användaren välja ett område att zooma in i genom att markera det önskade layoutområdet. Klicka på höger musknapp för att zooma ut till 100 %.

## Plate Layout Window (Fönstret plattlayout)

I det centrala området av fönstret **Plate Layout (Plattlayout)** visas en schematisk layout av en mikroplatta.

När en ny metod skapas väljer du brunnarna för mätningen. Ett klick med musen väljer ut en individuell brunn. De omarkerade brunnarna förblir gråa.

För den definierade layouten visas de identifierare, transformationer och formler som valts i Kontrollfältet i motsvarande brunnar (se Kontrollfältet för Create/Edit Method (Skapa/redigera meny) nedan).

#### Plattlayout: Sammanhangsberoende meny

Genom att högerklicka på en brunn i plattlayouten visas en sammanhangsberoende meny för de markerade brunnarna. Följande kommandon är tillgängliga:

Meny	Beskrivning
Summary (Sammanfattning)	Any plate view. (Alla plattvyer) Dialogrutan Summary (Sammanfattning) visas. Se kapitel 7.4.18 Sammanhangsberoende meny för en brunn/Dialogrutan Summary (Sammanfattning) för ytterligare information om dialogrutan Sammanfattning. Det här alternativet är tillgängligt om en layout har tilldelats den utvalda brunnen.
Fill Selection (Fyll urval)	Endast i läget plattlayout: Om en brunn eller ett område på plattan har markerats kan det fyllas med respektive identifierare. ID- och färgidentifierare kommer att visas på plattlayouten.
Delete Selection (Radera urval)	Endast i läget plattlayout: Detta kommando raderar ID- och färgidentifierare för de utvalda brunnarna och lämnar dem blanka.
Select all unused (Välj alla oanvända)	Endast i läget plattlayout: Alla oanvända brunnar på plattan är markerade.
Set / Remove Alias… (Ställ in/ta bort alias)	<b>Endast i läget plattlayout</b> : Ställer in eller tar bort alias för brunnens namn. Se kapitel 4.3.2 Metodlayout: Hur man definierar en plattlayout/Tilldela alias till brunnar.

#### Kontrollfältet i fliken Create/Edit Method (Skapa/redigera metod)

I **Kontrollfältet** till vänster på skärmen finns ett antal alternativ som ska utföras enligt den föreslagna sekvensen. Beroende på typ av mätning och ansluten läsartyp kan det hända att vissa av alternativen inte är tillgängliga. De är därför dolda.

När du väljer ett alternativ i kontrollfältet visas motsvarande dialogruta eller plattvy i panelen **Plate Layout window (Fönstret plattlayout)** på höger sida av skärmen.

Alla tillgängliga alternativ markeras med ett kryss när de har definierats. För alla typer av transformationer visas namnet på transformationen.

Fliken Kontrollfält - Create/edit method (Skapa/redigera metod) innehåller följande element:

# 4. Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)

Gruppen <b>Method layout</b> (Metodlayout)	<ul> <li>Elementet Plate layout (Plattlayout)</li> <li>Fönstret Plate layout (Plattlayout) öppnas och visar plattlayouten och dialogrutan Well assignment (Tilldelning av brunn).</li> <li>Elementet Conc., Dil., and Refvalues (Koncentration, spädning och referensvärden)</li> <li>Dialogrutan Concentration/Dilution/Reference (Koncentration/spädning/referens) öppnas och visar spädning, koncentration eller referensvärde för varje brunn. Standardkoncentrationerna kan fastställas samt spädningsfaktorerna och referensvärdena. Funktionen Autofill (Autofyll) gör det enkelt att tilldela koncentrationer ifall det finns distinktiva matematiska samband mellan koncentrationerna i de individuella brunnarna.</li> </ul>	
Förhands- beräkning	Reducering av spektradata för våglängdsskanningar	
Gruppen <b>Transformed data</b> (Transformerad data)	<ul> <li>Elementet Add new transformation (Lägg till ny transformation)</li> <li>Fönstret Plate layout (Plattlayout) öppnas och visar plattlayouten och inmatningsfältet Transformations (Transformationer)</li> </ul>	
Gruppen <b>Kinetic</b> (Kinetik)	<ul> <li>Elementet Kinetic data reduction (Reducering av kinetiska data)</li> <li>Fönstret Kinetics Calculation Parameters (Kinetiska beräkningsparameterar) öppnas. Utvärderingen av data för kinetiska mätningar kan definieras.</li> </ul>	
Kinetics transformation (Kinetiska transformationer)	• Elementet Add new kinetic transformation (Lägg till ny kinetisk transformation) Fönstret Plate layout (Plattlayout) öppnas och visar plattlayout och inmatningsfältet Transformations (Transformationer) (endast kinetiska indata kan väljas).	
Gruppen <b>Concentrations</b> (Koncentrationer)	• Elementet <b>Standard curve (</b> Standardkurva) Fönstret <b>Standard Curve</b> (Standardkurva) öppnas för att ställa in parametrarna för beräkning av koncentrationer och diagramvisning av standardkurvan för utvärderade data.	
Gruppen Concentration transformation data Koncentrations- transformationer)	<ul> <li>Elementet Add new concentration transformation (Lägg till ny koncentrationstransformation)</li> <li>Fönstret Plate layout (Plattlayout) öppnas och visar plattlayouten och inmatningsfältet Concentration Transformations (Koncentrationstransformationer) (endast koncentrationer kan väljas som indata).</li> </ul>	
Gruppen <b>Evaluate data</b> (Utvärdera data)	<ul> <li>Elementet Cutoff definition (Definiton av cutoff)</li> <li>Elementet QC validation (QC-validering) Fönstret Define Cutoff (Definiera cutoff) öppnas. I den här dialogrutan kan gränserna för en kvalitativ utvärdering (screening) definieras.</li> <li>Fönstret Define QC Validation (Definiera QC- validering) öppnas. Det används för att verifiera giltigheten hos ett test.</li> </ul>	

Gruppen <b>Data handling</b> (Datahantering)	<ul> <li>Elementet Data export (Dataexport)</li> <li>Elementet Printed report (Utskriven rapport)</li> <li>Elementet Automated data handling (Automatiserad datahantering)</li> </ul>
Gruppen <b>Miscellaneous</b> (Övrigt)	<ul> <li>Elementet User prompts (Användarprompt)</li> <li>Elementet Number format (Talformat)</li> <li>Elementet Method notes (Metodanteckningar)</li> <li>Fönstret <b>Define User Prompts</b> (Definiera användarprompt) gör det möjligt att tilldela data (nyckelord, kommentarer eller prompter) för varje mätning, som sedan kan tas med i en utskrift.</li> </ul>
	I fönstret <b>Number format</b> (Talformat) kan användaren definiera talformatet för rådata eller transformerade data som visas. I <b>Method notes</b> (Metodanteckningar) är det möjligt att ange en beskrivning av metoden.

# Kontrollfältet – fliken Create/Edit Method (Skapa/redigera metod) Sammanhangsberoende meny

Genom att högerklicka på en transformation, kinetisk transformation och koncentrationstransformation i kontrollfältet visas en sammanhangsberoende meny för den markerade transformationen. Följande kommandon är tillgängliga:

Meny	Beskrivning
Rename transformation (Byt namn på transformation)	Ett annat namn kan tilldelas den valda transformationen.
Insert transformation (Infoga transformation)	Används för att definiera en ny transformation.
Remove transformation (Ta bort transformation)	Används för att ta bort en transformation.

# 4.3.2 Metodlayout: Hur man definierar en plattlayout

Expandera Method layout (Metodlayout) i Kontrollfältet och välj Plate layout (Plattlayout). Plate view (Plattvyn) och dialogrutan Well Assignment (Tilldelning av brunn) visas.

Varje brunn i analysplattan kan tilldelas en identifierare. På grund av detta bör plattlayouten definieras (dvs. en brunn som anses vara positiv kontroll ska tilldelas en annan identifierare än en brunn som anses vara negativ kontroll).



Standardidentifierare är:

/ell Assignm.		Sample (Prov)	SM (Sample - Prov)
Identifiers SM BL	Exp. group:	Blank	<b>BL</b> (Blank) <b>BF</b> (Polarization reference buffer - Polarisering för referensbuffer)
ST PC		Reference (Referens)	<b>RF</b> (Polarization reference - Polariseringsreferens)
NC		Standard	ST (Standard)
LPC	Replicates	Control (Kontroll)	PC (Positive control - Positiv kontroll)
BF			<b>NC</b> (Negative control - Negativ kontroll)
RF			<b>LPC</b> (Low positive control - Svagt positiv kontroll)
Def. identif.			<b>HPC</b> (High positive control - Starkt positiv kontroll)
Fill selection	Delete selection		CL (Calibrator - Kalibrator)

## Tilldela en identifierare till brunnen

I dialogrutan **Well Assignment (Tilldelning av brunn)** kan identifiering av brunnar och layoutdefinition utföras. Den här lådan har också ett antal automatiserade funktioner för tilldelning av ID, vilket är ett viktigt verktyg för plattor av hög densitet.

De önskade inställningarna måste göras i dialogrutan **Well Assignment** (Tilldelning av brunn). Vanligtvis kan brunnarna markeras genom att klicka på en individuell brunn eller dra musen över de önskade brunnarna.

Följande sätt är möjliga för att tilldela brunnarna den valda identifieraren:

- Dubbelklicka när du gör urvalet av brunnen.
- Välj brunnarna på mikroplattan och klicka sedan på knappen Fill selection (Fyll urval) (eller klicka på höger musknapp och välj Fill selection (Fyll urval) i den sammanhangsberoende menyn) i dialogrutan Well Assignment (Tilldelning av brunn).
- Välj brunnarna på mikroplattan och dubbelklicka sedan på identifieraren i listan för identifierare i dialogrutan för tilldelning av brunn.
- När du har tilldelat definitionerna ändras displayen i de ändrade brunnarna.

#### Exempel på en märkt brunn:

SM1_4	1:a raden: prov, försöksgrupp nummer 1, prov-ID nummer 4.
1/14	2:a raden: antal replikat är 1, det totala antalet replikat är 14.
x-BL1	3:e raden: primärt tom - fylls med transformationsformeln, eller ett koncentrations-, spädnings-, referensvärde, om det är definierat, t.ex. minska värdet för brunnen (x) med medelvärdet av blanka värden.

Grupprutan Identifiers (Identifierare)	<ul> <li>I denna gruppruta ska motsvarande identifierare för de markerade brunnarna väljas.</li> <li>Alla tillgängliga identifierare visas i en rullgardinsmeny.</li> <li>Klicka på knappen <b>Define Identif</b> (Definiera identifierare) för att definiera ytterligare identifierare. Dialogrutan <b>Define Identifiers</b> (Definiera identifierare) visas, se nedan för mer information.</li> </ul>
Urvalsfältet Exp. group (Försöksgrupp)	Om plattan består av mer än ett test måste det finnas mer än en försöksgrupp. I urvalsfältet <b>Experiment group</b> (Försöksgrupp) definierar du vilket experiment brunnarna tillhör.
Grupprutan ID-Num. (ID-nummer)	<ul> <li>ID Number (ID-nummer) används för att tilldela samma ID till replikat som hör samman. ID-numret är endast tillgängligt för prover och standarder.</li> <li>Urvalsfältet ID-Num. (ID-nummer) Alternativet ID-Num (ID-nummer) tillåter användaren att markera de replikat som är associerade med samma ID. Detta kan användas om replikaten har tilldelats olika områden på plattan.</li> <li>Pilknappar ID-numret beräknas automatiskt. Om ett antal brunnar är markerade kan man med hjälp av pilknapparna bestämma i vilken riktning ID:n tilldelas brunnarna (vertikalt, horisontellt).</li> </ul>
Grupprutan Replicates (Replikat)	<ul> <li>Bestämmer antalet replikat för den valda identifieringstypen. Två alternativknappar gör det möjligt att bestämma om flera eller individuella värden ska definieras:</li> <li>Alternativknappen Fix number (Fast nummer) Aktiveras endast för standarder och prover där ID:n kan användas. Om denna alternativknapp är aktiv kan ett nummer anges i motsvarande textfält. Detta nummer anger hur många replikat som är avsedda för denna identifierare. De utvalda brunnarna fylls sedan med det angivna antalet replikat. Därför måste antalet valda brunnar vara en multipel av det angivna antalet replikat.</li> <li>Alternativknappen AII (Alla) Alla utvalda brunnar definieras som replikat för en identifierare. Om ett befintligt ID-nummer för proverna och standarderna väljs, läggs de utvalda brunnarna till som replikat till de befintliga replikaten. Med alla andra identifieringstyper läggs de utvalda brunnarna till som replikat till de befintliga replikaten.</li> </ul>
Knappen Fill section (Fyll urval)	Om ett område på plattan har markerats kan det fyllas med respektive identifierare. ID:n och färgen på identifierarna kommer att visas på plattlayouten.
Knappen Delete section (Radera urval)	Klicka på <b>Delete (Radera urval)</b> eller tryck på <b>DEL</b> för att radera ID:n och färgidentifierare för de utvalda brunnarna, och lämna dem blanka.

Dialogrutan Well assignment (Tilldelning av brunn) innehåller följande element:



KnappenKlicka på Def. identif (Definiera identifierare) om en ny<br/>identifierare ska definieras eller om en befintlig identifierare<br/>ska redigeras (se Expertens kunskap, Definiera en ny<br/>identifierare).

#### Tilldela alias till brunnar

Om du vill tilldela alias för definierade namn på brunnar klickar du på höger musknapp i den önskade brunnen, välj **Set/Remove Alias... (Ställ in/ta bort alias)** och välj ett identifieringsnamn i rullgardinsmenyn. Aliasnamnet är markerat med en asterisk \* och tillhör samma försöksgrupp, har samma ID-nummer och replikatnummer som den primärt definierade brunnen. Denna funktion används om t.ex. 0-standarden också används som negativ kontroll.

#### Definiera en ny identifierare

l dialogrutan Well Assignment (Tilldelning av brunn), klicka på Def. Identif för att definiera en ny identifierare eller redigera en befintlig.

Constant and the second	Identifier		
Identifier	Type:	Sample	
SM	Abbreviati	on: SM	Cancel
Blank BL BF BF BF BF BF BF BF Standard ST	Descriptio	n: Sample	Help
Control PC NC LPC	<b>~</b>	ave Discard	Set as default
Edit			Jet as derault
Delete			Bestore defaul

Identifierarna är listade i grupper (se tabellen nedan). Om du markerar en identifierare visas dess egenskaper i det högra fönstret.

Standardidentifierare är:

Sample (Prov)	SM (Sample - Prov)
Blank	BL (Blank)
	<b>BF</b> (Polarization reference buffer - Polarisering för referensbuffer)
Reference (Referens)	<b>RF</b> (Polarization reference - Polariseringsreferens)
Standard	ST (Standard)
Control (Kontroll)	<ul> <li>PC (Positive control - Positiv kontroll)</li> <li>NC (Negative control - Negativ kontroll)</li> <li>LPC (Low positive control - Svagt positiv kontroll)</li> <li>HPC (High positive control - Stark positiv kontroll)</li> <li>CL (Calibrator - Kalibrator)</li> </ul>



Trädstrukturen för Identifier (Identifierare)	En strukturerad vy av alla befintliga identifierare, deras färger och förkortningar visas i ett litet fönster. Identifierarna är listade under grupperna <b>Sample</b> (Prov), <b>Blank, Reference</b> (Referens), <b>Standard</b> (Standard) och <b>Control</b> (Kontroll).
Grupprutan Identifier (Identifierare)	Kriterierna för de olika identifierare som används i programmet visas. Om nya identifierare behövs kan de läggas till här.
	I grupprutan <b>Identifier (Identifierare)</b> visas typ, förkortning, färg på brunn och beskrivning av den identifierare som valts i trädstrukturen.
	Grupprutan <b>Identifier (Identifierare)</b> aktiveras endast för att ange data när
	<ul> <li>en av identifierargrupperna i trädstrukturen är markerad och knappen Insert (Infoga) har klickats eller när</li> </ul>
	<ul> <li>en av identifierarna i trädstrukturen är markerad och knappen Edit (Redigera) har klickats.</li> </ul>
	Därefter kan du i den här grupprutan definiera typ, förkortning, färg på brunn och beskrivning av den nya identifieraren.
	Textfältet <b>Type (Typ)</b> : Typ av identifierare visas. Inga ändringar är möjliga.
	Textfältet <b>Abbreviation (Förkortning)</b> : Identifieringens förkortning kommer att användas i visningen av analysplattan.
	Rullgardinsmenyn <b>Color (Färg)</b> : Färgen på identifieraren på plattan måste väljas här.
	Textfältet <b>Description (Beskrivning)</b> : En textbeskrivning kan anges för varje identifierare.
	Använd en av dessa två knappar för att spara ändringar av nyligen angiven data eller för att avvisa eventuella ändringar:
	Knappen <b>Save (Spara)</b> : Knappen <b>Save (Spara)</b> sparar den angivna färgen, förkortningen och beskrivningen av den relevanta identifieraren.
	Knappen <b>Discard (Avvisa):</b> Knappen Discard (Avvisa) avbryter alla ändringar.
Knappen <b>Insert (Infoga)</b>	Klicka på knappen <b>Insert (Infoga)</b> för att skapa en ny identifierare . Denna nya identifierare kommer att associeras med den aktuellt valda identifierargruppen i trädstrukturen.
Knappen <b>Edit (Redigera)</b>	En vald identifierare kan ändras.
Knappen <b>Delete (Radera)</b>	Klicka på <b>Delete (Radera)</b> eller tryck på <b>DEL</b> för att radera alla utvalda identifierare.

Dialogrutan **Define Identifiers (Definiera identifierare)** innehåller följande element:



Knappen Set as default (Ställ in som standard)	Med det här alternativet kan inställningarna definieras som standardinställningar för framtida användning.
Knappen Restore default (Återställ till standard)	Med det här alternativet kan inställningarna återställas till de tidigare definierade standardvärdena.

# 4.3.3 Metodlayout: Koncentration, spädning och referensvärden

I Kontrollfältet, expandera Method layout (Metodlayout) och välj Conc./Dil./Ref.-values (Koncentration/spädning/referensvärden). Plate view (Plattvyn) och dialogrutan Select Identifier (Välj identifierare) visas.



#### Dialogrutan Concentration/Dilution/Reference

(Koncentration/spädning/referens) innehåller följande element:

Rullgardinsmenyn <b>Select Identifier</b> (Välj identifierare)	Värdena är relaterade inom de individuella brunnstyperna. Detta fält anger alla brunnar som för närvarande ingår i mätningen. Välj en identifierare för att redigera motsvarande spädnings- eller koncentrationsfaktorer. Alla brunnar som matchar den utvalda brunnstypsidentifieraren listas.
Urvalsfältet <b>Exp. group</b> (Försöksgrupp)	Välj respektive försöksgrupp. Om det bara finns en försöksgrupp kan fältet inte redigeras.

Tabell med kolumnerna <b>Identifier</b> (Identifierare) och <b>Concentration/</b> <b>Dilution</b> (Koncentration/ spädning).	Tabellen visar alla brunnar i kolumnen <b>Identifier</b> (Identifierare) som matchar den utvalda identifieraren (från listan <b>Select Identifier</b> (Välj identifierare)). Motsvarande spädningar, koncentrationer eller referensvärden infogas och listas i kolumnen <b>Dilution/Concentration/Reference</b> (Koncentration/spädning/referens). Om du t.ex. anger en spädningsfaktor på 2 betyder det att provet har spätts ut med hälften. Den beräknade koncentrationen multipliceras därför med 2.
Textfältet Unit (Enhet)	Den visade koncentrationsenheten kan bestämmas.
Kryssrutan <b>Autofill</b> (Autofyll)	Funktionen <b>Autofill</b> (Autofyll) gör det möjligt att beräkna motsvarande koncentrationer eller spädningsfaktorer enligt tillgängliga typer av serier.
Knapparna <b>Option</b> (Alternativ) för matematisk koncentrations- beräkning	<ul> <li>Om kryssrutan Autofill (Autofyll) har markerats finns följande alternativ tillgängliga:</li> <li>Aritmetisk talföljd: a (n+1) = a (n) +</li> <li>Geometrisk talföljd: a (n+1) = a (n) x</li> <li>Användardefinierad talföljd a (n+1) = Exempel: Varje efterföljande koncentration ska innehålla två gånger + 0.5 av den föregående koncentrationen: Formel 1: a(n+1) = 2 * n + 0.5</li> </ul>
Knappen <b>Apply</b> (Tillämpa)	Knappen <b>Apply</b> (Tillämpa) tillämpar den valda matematiska koncentrationsberäkningen på brunnarna som visas i tabellen med kolumnerna <b>Identifier</b> (Identifierare) och <b>Concentration/ Dilution</b> (Koncentration/spädning)
Grupprutan <b>Dilution series</b> (Spädningsserie)	Kryssrutan <b>Calculate ICx</b> (Beräkna ICx) Visar ett spädningsdiagram för provet och beräknar automatiskt ICx-värdena. Detta kräver prover med minst 4 replikat och minst 4 olika spädningar.
	Rullgardinsmenyn <b>Input Data</b> (Indata) Välj indata från rullgardinsmenyn.
	Textfältet <b>Calculation Condition</b> (Beräkningsvillkor) Skärningspunkten beräknas med den angivna procentandelen av det maximala värdet respektive det maximala värdet minus det minsta värdet*.
	Textfältet <b>ICx name</b> (ICx-namn) Data kan väljas för beräkning av ICx. Namnet fylls i automatiskt baserat på beräkningsvillkoret.
	<b>Set 0% value to</b> (Ställ in 0 %-värdet till) Intensitet 0* Min. intensitet i spädningsserien*

Spädningen kan redigeras direkt i plattlayouten genom att högerklicka på en eller flera utvalda brunnar. På detta sätt är det möjligt att tilldela replikaten olika spädningsvärden.

## Beräkning av ICx

Den matematiska beräkningen av anpassningen för spädningsserien är identisk med beräkningen av standardkurvan med Marquardts algoritm med fyra parametrar.

Detta kräver minst fyra replikat med olika spädningar.

Dessutom beräknas det/de specificerade skärningspunkten/skärningspunkterna t.ex. IC50.



Det är möjligt att definiera fler än en skärningspunkt för spädningsserien.

Det är möjligt att definiera 0 %-värdet för beräkningen genom att välja:

Intensitet 0

Använda 0 OD

Det största värdet i spädningsserien anses vara 100 %, värdet 0 anses vara 0 %. ICx (t.ex. IC50) definieras sedan som den spädning där responsen når x % (t.ex. 50 %). Det beräknas endast om värdet ligger inom tillgängliga data (ingen extrapolering används).

eller

Lägsta intensitet i spädningsserien
 Användning av den lägsta intensiteten i spädningsserien



#### Anmärkning

IC50 beräknas ofta med medelvärdet för spädning = 1:1 och spädning = 1:oändligt. För att uppnå detta måste spädningsserien vara korrigerad för utgångsvärdet före beräkningen av ICx.

Det största värdet i spädningsserien anses vara 100 %, den minsta intensiteten i spädningsserien anses vara 0 %.



# 4.3.4 Förhandsberäkning: Reducering av spektradata

Det här alternativet är endast tillgängligt för mätningar som innehåller en tvådimensionell skanning (absorbansskanning).

Expandera **Precalculation (Förhandsberäkning)** i kontrollfältet och välj **Spectra Data Reduction (Reducering av spektradata)**.

Reducering av spektradata gör det möjligt att utföra en spektrablankreducering och extrahera specifika data, t.ex. intensiteter från skanningen.

# Dialogrutan **Spectra data reduction (Reducering av spektradata)** innehåller följande element:

Data reduction		Available results:
Blank reduction	Options	
Smooth	Options	S Smoothed spectra
🗹 Wavelength pick	Options	S Normalized spectrum
🗹 Normalize	Options	V Peak intensity V Wavelength peak
🗹 Peak	Options	V Peak width
Custom formulas	Options	Y Peak area
Aryssrutan Hank reduction Blank- educering)	har defin beräknas från alla på plattar ytterligar definiera	ierats i layouten. Spektrablankreduceringen s genom att subtrahera blankbrunnens spektrum andra brunnar. När fler blanka värden definieras n, t.ex. när flera försöksgrupper definieras, visas e en alternativknapp. Detta gör det möjligt att det blanka värdet som ska användas.
	Textrutar behandla	n <b>Input data (Indata)</b> : visar de indata som ska Is.
	Alternativ välj om e försöksgi	/knappen <b>Blank reduction (Blankreducering)</b> : n blankreducering ska utföras på den utvalda ruppen.
	Rullgardi välj för at för blank	nsmenyn <b>Reduce all by (Reducera alla med)</b> : It definiera den identifierare som ska användas reducering.
Kryssrutan Smooth Utjämning)	Ett spekt En utjäm faktor de	rum som innehåller mycket brus kan jämnas ut. ningsfaktor kan ställas in bland alternativ. Denna finierar graden av utjämning.
	Textrutar behandla	n <b>Input data (Indata)</b> : visar de indata som ska is.
	Smooth utjämning	factor (Utjämningsfaktor): för att definiera gsfaktorn.



Kryssruta för val av våglängd	Wavelength pick (Val av våglängd) används för att extrahera intensiteter vid specifika våglängder och för beräkna resultat som kvoter. Det är obligatoriskt att stä in alternativen.	att illa		
	Dialogrutan <b>Wavelength pick (Val av våglängd)</b> innehåller följande element:			
	Wavelength pick parameters	Wavelength pick parameters X		
	Input data: Smoothed spectra			
	○ Pick Wavelengths       WL [nm]       500       ✓         ○ Calculate Ratio       WL [nm]       600       ✓       /       WL [nm]       700       ✓         ○ Area       from WL [nm]       500       ✓       to WL [nm]       700       ✓         ○ Custom       Formula			
	Wavelengths     Constants     Operators       500       +			
	Data name         Formula         Add           500 nm         x(500)         Delet           600 nm / 700 nm         x(600)/x(700)         Edit	e		
	< >> OK Cancel Help	)		
	<ul> <li>anges.</li> <li>Textrutan Input data (Indata): visar de indata som ska behandlas.</li> <li>Alternativknappen Pick Wavelengths (Välj våglängde en intensitet vid en specifik våglängd kan läggas till i lis av beräknade resultat.</li> </ul>	a er): stai		
	Alternativknappen <b>Calculate Ratio (Beräkna kvot)</b> : tv våglängder kan anges. Kvoten mellan intensiteterna vi dessa våglängder beräknas och finns tillgängliga som resultat.	å d		
	Area: för att beräkna arean under spektrakurvan mella två definierade våglängder.	เท		
	Alternativknappen <b>Custom (Anpassad)</b> : en användardefinierad formel kan anges. Intensiteter vid specifika våglängder kan användas i formler.			
	Listrutan <b>Results (Resultat)</b> : sammanfattar alla definierade formler.			
	Knappen <b>Add (Lägg till)</b> : det aktuella urvalet läggs till resultatlistan.	i		
	Knappen <b>Delete (Radera)</b> : det valda resultatet tas bor från listan.	t		
	Knappen <b>Edit (Redigera)</b> : datanamnet för det valda resultatet kan definieras.			

# 4. Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)

Kryssrutan Normalize	Dialogrutan Normalize parameters (Normalisera parametrar) innehåller följande element:				
(Normalisera)	Normalize parameters X				
	Input data: Smoothed spectra				
	Normalize to Cancel				
	Max. intensity				
	O Intensity at WL [nm] 500 Help				
	Använd det här alternativet för att normalisera spektrumet till intensiteten vid en viss våglängd eller till den maximala intensiteten i spektradiagrammet för varje brunn.				
	Textrutan <b>Input data (Indata)</b> : visar de indata som ska behandlas.				
	Normalize to (Normalisera till): välj mellan maximal intensitet eller infoga anpassad intensitet.				
Kryssrutan Peak (Topp)	En topp definieras som positionen för det maximala värdet i spektrumet. Överflöden och maxima vid spektrumets vänstra eller högra kant behandlas inte som toppar. Det är mycket viktigt att spektrumet är tillräckligt jämnt innan man letar efter toppar.				
	Dialogrutan <b>Peak parameters (Parametrar för toppar)</b> innehåller följande element:				
	Peak parameters X				
	Input data: Normalized spectrum OK				
	Wavelength range: from: 400 nm Cancel				
	to: 600 nm				
	Ignore peaks below 0 OD/RFU				
	Calculate width and area at 50 % of max. intensity Help				
	Använd det här alternativet för att hitta toppen med den högsta intensiteten inom en definierad våglängd.				
	Tröskelvärdet (OD/RFU-värde) och beräkningskriterierna för beräkning av bredd och area kan också definieras.				
	Följande data beräknas för den funna sökvägen:				
	toppens intensitet				

- toppens våglängd
- toppens bredd
- toppens område



Kryssrutan Custom formulas (Anpassade formler)	Markera den här kryssrutan om du vill ange formler för att beräkna spektrumet med hjälp av givna funktioner, t.ex. utjämning. Se kapitel 11.2 Reducering av spektradata för ytterligare information.
	Transformation måste antingen ge ett spektrum eller en enda data för alla brunnar; det är inte möjligt att kombinera båda. Med det här alternativet kan man definiera funktioner som inte är tillgängliga i standardurvalet. T.ex. för att hitta fler toppar i ett spektrum eller för att skapa en härledning av ett spektrum.
Tillgängliga resultat	Tillgängliga resultat för alla data listas. S är ett spektrum V är ett värde.

# 4.3.5 Transformed data (Transformerad data): Lägg till ny transformation

l kontrollfältet, expandera **Transformed data (Transformerad data)**. Alla definierade transformationer visas i kontrollfältet.

Om du vill definiera en ny transformation väljer du den eller de brunnar som transformationen ska tillämpas på och klickar på Add new transformation... (Lägg till ny transformation).

Standardnamnet bör ändras omedelbart i redigeringsrutan. Namnet kan infogas eller ändras senare genom att klicka på **Rename Transformation (Byt namn på transformation)** i den sammanhangsberoende menyn.

Ett typiskt exempel på en transformation är reduceringen av det tomma värdet (blank) från alla brunnar.

#### Anmärkning

Namnet på transformationen måste vara i latinska bokstäver för att göra transformationen tillgänglig som indata för ytterligare transformationer med flera indata.

#### Anmärkning

Namnet på en transformation används för att representera resultatet av beräkningarna och kommer också att användas när värdena visas i de specifika menyerna för datautmatning. Beräknade värden för transformationerna finns också tillgängliga som indata för ytterligare utvärderingar.



#### Anmärkning

Om ett blankt värde har ställts in i plattlayouten föreslås formeln för att beräkna blankreduceringen i kombinationsrutan: x-BL1.

Symbolen x hänvisar till det aktuella värdet i en brunn. BL1 är medelvärdet för blankbrunnen/blankbrunnarna.



#### Anmärkning

Om en mätning med multipla inmärkningar med två märkningar definieras, föreslås olika formler för att reducera data med multipla inmärkningar i kombinationsrutan.

'Label1'!x/'Label2'!x ... fördefinierad kvotberäkning 'Label2'!x/'Label1'!x ... fördefinierad kvotberäkning 'Label1'!x-'Label2'!x ... fördefinierad differensberäkning ('Label2'!x-'Label1'!x ... fördefinierad differensberäkning ('Lable1'!x-'Label1'!BL1)/('Label2'!x-'Label2'!BL1) ... fördefinierad kvotberäkning med blankreducering ('Lable2'!x-'Label2'!BL1)/('Label1'!x-'Label1'!BL1) ... fördefinierad kvotberäkning med blankreducering



#### Anmärkning

Om läsningsläget är inställt på absorbans föreslås formeln för att beräkna transmissionen i kombinationsrutan: 1/10<sup>x</sup>.

Ett antal transformationsberäkningar kan definieras när metoden konfigureras. Ett typiskt exempel på detta är borttagning av det tomma värdet (blank) från alla brunnar. Se kapitel 11.3 Hur man skriver en formel för mer detaljer.

Textrutan för formler ovanför plattvyn används för att definiera olika transformationsberäkningar, med hjälp av rådata och tidigare transformationer som indata.

På mikroplattan måste man välja de brunnar där transformationen ska tillämpas.

Redigeringsfältet för transformationer innehåller följande element:

Rullgardinsmenyn Input Data (Indata)	Listan innehåller mätvärdena, resultaten från förhandsberäkningar och alla redan definierade transformationer samt medelvärden. Om data för <b>Mean</b> (Medelvärde) väljs som Input data (Indata) för Transformations (Transformationer) kommer transformationerna endast att tilldelas det första replikatet av en identifierare.
Textrutan Formula (Formel) fx	Formler kan anges genom att skriva in eller genom att välja den funktion som behövs från listrutan Functions&Constants (Funktioner och konstanter).
Rullgardinsmenyn <b>Formula (Formel)</b>	Listan innehåller vissa standardformler (se Anmärkningar nedan) och alla formler som har inkorporerats i den nuvarande metoden. En formel kan därför väljas från denna lista eller så kan en ny formel kan läggas till. Formler för transformationer kan anges med hjälp av lämpliga variabler, operatorer och multipla funktioner.
Knappen <b>Confirm (Bekräfta)</b>	Endast i redigeringsläget för formler. Grön bock Tilldelar transformationerna till den utvalda brunnen. Klicka på knappen <b>Confirm (Bekräfta)</b> eller tryck på <b>ENTER</b> för att tilldela formeldefinitionen till brunnen och ändra till läget <b>Select (Välj)</b> .



Knappen <b>Cancel (Avbryt)</b>	Endast i redigeringsläget för formler. Rött kryss Klicka på knappen <b>Confirm (Bekräfta)</b> eller tryck på tangenten <b>CANCEL</b> för att lämna läget <b>Edit (Redigera)</b> utan att tilldela formeldefinitionen till brunnen.
Rullgardinsmenyn Available data (Tillgängliga data)	Välj lämpliga data från listan om mer än ett set indata ska användas för beräkningar. Datasetet visas i formelns textruta inom apostrofer följt av ett utropstecken. För att slutföra definitionen anger du identifierarens namn eller hänvisar till motsvarande värde i brunnen. T.ex. 'Rådata' ! BL1 Listan innehåller mätvärden, resultat från förberäkningar och alla hittills definierade transformationer.
Rullgardinsmenyn Functions (Funktioner)	Listar alla matematiska och booleska funktioner för att definiera formler.
Knappen <b>Options (Alternativ)</b>	<ul> <li>Använd den här knappen för att anpassa beteendet när du väljer brunnar i redigeringsläget. Det går att välja mellan följande alternativ:</li> <li>Identifierare genomsnittsnamn</li> <li>Identifierare replikatnamn</li> <li>Namn på plattans brunn</li> </ul>
Knappen Constants (Konstanter)	Klicka på den här knappen för att öppna dialogrutan <b>Define Constants (Definiera konstanter)</b> där konstanter för beräkningar kan definieras.

Definitionen för transformation har två lägen:

Select mode (Urvalsläge)	När du klickar på en brunn visas motsvarande formel i textrutan. Den aktiva brunnen visas med en röd kant.
Edit mode (Redigeringsläge)	När du skriver in en formel eller trycker på tangenten "=" visas den aktiva brunnen med en blå kant. Om du väljer andra brunnar läggs motsvarande identifierare till den aktuella formeln. När du väljer den aktiva brunnen läggs ett "x" till i formeln. Symbolen "x" avser det aktuella värdet i en brunn.

När formeln har skapats tilldelas den till den utvalda brunnen om du trycker på enter eller klickar på bekräftelseknappen (grön bock). Efter att ha tilldelat formeln nås urvalsläget. En integrerad formelkontrollant verifierar att formeln är trovärdig. En varning visas om den skapade formeln inte är möjlig att genomföra.

Välj **CANCEL (Avbryt)** för att avvisa den definierade formeln och nå det valda läget. Det är också möjligt att använda motsvarande knappar till vänster om formeltextrutan.

I läget **Select mode (Urvalsläge)** klickar du på den röda rutan i brunnens högra nedre kant och drar den röda urvalsramen över de brunnar där formeln ska gälla.

Det är också möjligt att välja önskade brunnar och använda funktionerna **Set Formula (Ställ in formel)** och **Remove Formula(s) (Ta bort formel/formler)** som är tillgängliga från den sammanhangsberoende menyn.

Dessa åtgärder måste utföras för alla brunnar som ska innehålla transformationer.



#### Definiera konstanter

Använd dialogrutan **Define Constants (Definiera konstanter)** för att definiera konstanta värden i en metod. Dessa konstanter kan användas överallt där en formel kan anges

Klicka på knappen **OK** för att spara de ändrade parametrarna.

Dialogrutan **Define Constants (Definiera konstanter)** innehåller följande element:

Listan <b>Constants</b> (Konstanter)	I varje rad kan en konstant definieras. Raderna är indelade i fyra kolumner:
	<ul> <li>Name (Namn) Ange ett lämpligt identifieringsnamn för konstanten, t.ex. en kod eller en förkortning. Endast bokstäver får användas.</li> </ul>
	<ul> <li>Value (Värde) I textfältet Value (Värde) måste ett numeriskt värde tilldelas konstanten.</li> </ul>
	Comment (Kommentar)     Skriv en kort kommentar om konstanten.
	<ul> <li>Req. (Krav) ("required") Kryssrutan Req. (Krav) anger att en mätning endast kan påbörjas när värdet för konstanten har mottagits eller bekräftats.</li> </ul>

## 4.3.6 Kinetik: Reducering av kinetiska data

l Kontrollfältet, expandera Kinetic (Kinetik) och välj Kinetic data reduction (Reducering av kinetiska data).

Dialogrutan är strukturerad i flikar: **Slopes**, **Onsets**, **Min./Max./Area**, **Available output data**, **Enzyme kinetics**. Enheten för värdena visas i enlighet med det valda mätläget (t.ex. OD för absorbans).



Anmärkning Oanvändbara data (t.ex. överflödsvärden) ignoreras vid beräkning av kinetiska data.

## Fliken Slopes (Lutning)

Den här fliken tillåter användaren fastställa utvärderingen av den kinetiska kurvans lutning:

Rullgardinsmenyn Input Data (Indata)	Välj indata som ska behandlas.
Rullgardinsmenyn <b>Calculation</b> Beräkning)	Välj beräkningsmetod, linjär eller kvadratisk. (se kapitel 11 Beräkningar).
Kryssrutan <b>Mean slope</b> (Genomsnittlig lutning)	Start- och sluttid kan skrivas in i de angivna textfälten, annars analyseras hela kinetiken.

	<ul> <li>Alternativknappen Time/Points (Tid/punkter) Om Time (Tid) är valt, anges start- och sluttiden i timmar, minuter och sekunder. Om Points (Punkter) är valt anges start- och slutpunkten för analysen genom det angivna numret för kinetikcykeln.</li> </ul>
	<ul> <li>Fältet Start</li> <li>Starttiden (i timmar, minuter och sekunder) eller cykelnumret måste anges här.</li> </ul>
	<ul> <li>Fältet End (Slut) Sluttiden (i timmar, minuter och sekunder) eller cykelnumret måste anges här.</li> </ul>
Kryssrutan Maximum slope (Maximal lutning)	Start- och sluttid kan skrivas in i de angivna textfälten, annars analyseras hela kinetiken.
	<ul> <li>Alternativknappen Time/Points (Tid/punkter) Om Time (Tid) är valt, anges start- och sluttiden i timmar, minuter och sekunder. Om Points (Punkter) är valt anges start- och slutpunkten för analysen genom det angivna numret för kinetikcykeln.</li> </ul>
	<ul> <li>Fältet Start Starttiden (i timmar, minuter och sekunder) eller cykelnumret ska anges här.</li> </ul>
	<ul> <li>Fältet End (Slut) Sluttiden (i timmar, minuter och sekunder) eller cykelnumret måste anges här.</li> </ul>
	<ul> <li>Textfältet <b>Points</b> (Punkter)</li> <li>Välj för hur många punkter beräkningen av maximal lutning ska utföras.</li> </ul>

Alternativet **mean slope (genomsnittlig lutning)** bestämmer den genomsnittliga lutningen (genomsnittlig ökning/minskning) för hela det definierade intervallet. En starttid, en sluttid och ett beräkningsläge (linjär eller kvadratisk) måste först anges. Alla mätpunkter som ligger inom det valda intervallet bestäms. En regressionslinje (linjär regression eller polynom av 2:a graden) läggs ut genom de valda mätpunkterna och den genomsnittliga lutningen bildas. Den genomsnittliga lutningen definieras som det aritmetiska medelvärdet av de lutningar som beräknas från mittpunkterna i två intilliggande mätpunkter. Den kinetiska data som görs tillgängliga med denna metod omfattar den genomsnittliga lutningen per sekund, minut och timme samt korrelationskoefficienten och anpassningsgraden.

Med alternativet **maximum slope (maximala lutningen)** bestäms den maximala lutningen (maximal ökning/minskning) som uppnås i det valda intervallet. Under alternativet **Points (Punkter)** ska antalet kombinerade punkter anges. I början beräknas lutningen från centrumpunkten för den 1:sta och den n:te punkten från de första n punkterna. Därefter flyttas intervallet en punkt ytterligare och processen upprepas. Denna process fortsätter över alla punkter inom det valda intervallet. Resultatet kommer att fastställas som det största absoluta värdet av dessa individuella lutningar. De kinetiska data som görs tillgängliga med denna metod omfattar den maximala lutningen per sekund, per minut och per timme samt tidsperioden från den första mätningen till den maximala lutningen i sekunder.

#### **Onsets (Starttider)**

Använd fliken **Onsets (Starttider)** för att bestämma tiden för att nå en viss datapunkt (OD-värde):

Rullgardinsmenyn Input Data (Indata)	Välj indata som ska behandlas.
Kryssrutan Time to onset (Tid till starttid)	Om kryssrutan <b>Time to onset</b> (Tid till starttid) är markerad kan du ange ett absolutvärde för starttidpunkten i följande textfält.
	<ul> <li>Textfält: Ett absolutvärde för starttiden måste anges.</li> </ul>
	<ul> <li>Alternativknapp och textfältet Basis mean of the first n points (Basmedelvärde av de första n punkterna): Ange önskat antal punkter om det är valt.</li> </ul>
	<ul> <li>Alternativknapp och textfält Basis: Om detta är valt måste ett absolutvärde för basen anges i det intilliggande textfältet.</li> </ul>
Kryssrutan <b>Time to onset %</b> (Tid till starttid %)	Om kryssrutan <b>Time to onset %</b> (Tid till starttid %) är markerad kan ett procentvärde för starttiden anges i följande textfält.
	<ul> <li>Textfält: Ett procentvärde för starttiden måste anges.</li> </ul>
	<ul> <li>Alternativknapp och textfältet Basis mean of the first n points (Basmedelvärde av de första n punkterna):</li> <li>Önskat antal punkter kan anges om det är valt.</li> </ul>
	<ul> <li>Alternativknapp och textfält Basis: Om detta är valt måste ett absolutvärde för basen anges i det intilliggande textfältet.</li> </ul>

De resultatdata som görs tillgängliga genom denna metod omfattar basvärdet, den tid det tar att nå basvärdet, den tid det tar att nå totalsumman av bas- och startvärdet och skillnaden mellan dessa två tidsperioder (Time Basis to Onset).

Med ökande kinetiska mätvärden ska startvärdet definieras som ett positivt tal, med minskande värden ska startvärdet definieras som ett negativt tal.

# Fliken Min./Max./Area

Använd den här fliken för att definiera utvärderingen av minsta och maximala värdena i kurvan och för att definiera beräkningen av arean under den kinetiska kurvan:

Rullgardinsmenyn <b>Input data</b> (Indata)	Välj indata som ska behandlas.
Kryssrutan <b>Mean minimum value</b> (Medelvärde av minsta värde)	Välj för att öppna textfältet <b>Points</b> (Punkter) där det fasta antalet punkter ska anges för det minsta värdet. En jämn kurva skapas sedan med hjälp av dessa punkter och det lägsta värdet i kurvan bestäms.



Kryssrutan <b>Mean maximum value</b> (Medelvärde av maximala värdet)	Välj för att öppna textfältet <b>Points</b> (Punkter) där det fasta antalet punkter ska anges för det maximala värdet. En jämn kurva skapas sedan med hjälp av dessa punkter och det högsta värdet i kurvan bestäms.
Kryssrutan <b>Area</b>	Start- och sluttid kan skrivas in i de angivna textfälten, annars analyseras hela kinetiken.
	<ul> <li>Alternativknappen Time/Points (Tid/punkter) Om Time (Tid) är valt, anges start- och sluttiden i timmar, minuter och sekunder. Om Points (Punkter) är valt anges start- och slutpunkten för analysen genom det angivna numret för kinetikcykeln.</li> </ul>
	<ul> <li>Fältet Start Starttiden (i timmar, minuter och sekunder) eller cykelnumret ska anges här.</li> </ul>
	<ul> <li>Fältet End (Slut) Sluttiden (i timmar, minuter och sekunder) eller cykelnumret måste anges här.</li> </ul>

Medelvärden tas från antalet punkter, med start från det första kinetiska mätvärdet. Detta pågår tills den sista kinetiska mätpunkten ingår i beräkningen för medelvärdet. Det minsta/maximala värdet bestäms utifrån dessa medelvärden.

De resultatdata som görs tillgängliga genom denna metod omfattar minsta/maximala värden och tidperioden från den första mätningen till det minsta/maximala värdet i sekunder.

#### Exempel:

Med en kinetik på 5 cykler och ett fast antal på 3 punkter tas följande medelvärden:

medelvärde av 1:a, 2:a och 3:e kinetiska värde

medelvärde av 2:a, 3:e och 4:e kinetiska värde

medelvärde av 3:e, 4:e och 5:e kinetiska värde

Det minsta/maximala värdet bestäms utifrån dessa tre medelvärden.

Arean under kurvan beräknas med hjälp av formeln nedan:

$$A = \sum_{i=1}^{n-1} y_i * (x_{i+1} - x_i) + \frac{(y_{i+1} - y_i) * (x_{i+1} - x_i)}{2}$$

# Fliken Available Data (Tillgängliga data)

En lista över resultaten visas på fliken Tillgängliga utdata:

Datafält

I datafältet visas de värden som valts i föregående flikar i dialogrutan Kinetic Calculation Parameters (Kinetiska beräkningsparameterar).

Detta datafält är endast avsett för visning av de angivna utdata och ger inga redigeringsfunktioner.



## Fliken Enzyme Kinetics (Enzymkinetik)

Använd den här fliken för att fastställa utvärderingen av enzymkinetik enligt modellen Michaelis-Menten:

Rullgardinsmenyn Input Data (Indata)	Välj indata som ska behandlas.
Kryssrutan <b>Calculate Km and Vmax</b> Beräkna Km och Vmax)	Bestäm om Km och Vmax ska beräknas genom att välja motsvarande kryssruta.
Grupprutan alculation type (Beräkningtyp)	Beräkningstypen kan väljas som:
	Hanes (koncentration mot koncentration/indata)
	Eadie-Hofstee (indata/koncentration mot indata)
	Lineweaver-Burk (1/indata mot 1/koncentration)

De resultatdata som görs tillgängliga genom denna metod omfattar Km och Vmax för enzymkinetiska diagram för varje försöksgrupp.

Till skillnad från resultaten av de andra kinetiska beräkningarna visas dessa resultat i dialogrutan Graph: Enzyme Kinetics (Diagram: Enzymkinetik).

Den här fliken är endast tillgänglig om standarder finns på plattan och om antingen transformationer eller lutningsberäkningar har definierats.

## 4.3.7 Kinetiska transformationer: Lägg till nya kinetiska transformationer

I dialogrutan Kinetic transformations (Kinetiska transformationer) kan en transformationsformel definieras. Formeln används för att transformera kinetiska indata individuellt för varje brunn.

I dialogrutan Kinetic transformations (Kinetiska transformationer) kan ytterligare beräkningar utföras på kinetiska indata.

Fönsterelementen är jämförbara med indata för transformation. Se kapitel 4.3.5 Transformed data (Transformerad data): Lägg till ny transformation för mer information.

# 4.3.8 Koncentrationer:Standardkurva

Använd det här alternativet för att ställa in standardkurvor för kvantitativa tester.

Dialogrutan för standardkurvan innehåller olika inställningar för analystyp, axlar och visning av standardkurvan. De redigerbara fälten och elementen är grupperade i 5 olika flikar.

# Fliken Data

Den här fliken innehåller några grundläggande inställningar, t.ex. källan för indata.

Spinnkontrollen	Om flera tester ska utföras på samma platta måste
Exp. group	Experiment group väljas. Om plattan endast innehåller
(Försöksgrupp)	ett test visas inte spinnkontrollen.
	För varje försöksgrupp kan indata definieras individuellt.



Rullgardinsmenyn Input Data (Indata)	Välj den <b>Input Data</b> (Indata) som ska användas för standardkurvan. Välj mätdata eller tillgängliga transformationsresultatet.
Alternativknappen Standards from Iayout (Standarder från layouten)	Beräkna standardkurvan från standarderna på layouten.
Alternativknappen Standards from ext. file (Standarder från extfilen	Om den valda försöksgruppen inte innehåller några standarder kan en standardkurva laddas från en .std-fil. Du måste klicka på knappen <b>Select</b> (Välj) för att välja filen.
Alternativknappen Standards from exp. group (Standarder från försöksgruppen)	Om den valda försöksgruppen inte innehåller några standarder kan en standardkurva från en annan försöksgrupp användas.
Alternativknappen <b>No standard curve</b> (Ingen standardkurva)	Om den valda försöksgruppen inte innehåller några standarder kan man besluta att inte beräkna några koncentrationer för denna försöksgrupp (i detta fall inställd som standard).
Knappen Additional Concentrations (Ytterligare koncentrationer)	Klicka på knappen <b>Additional Concentrations</b> (Ytterligare koncentrationer) för att öppna dialogrutan <b>Calculate Additional Concentration</b> (Beräkna ytterligare koncentrationer). Välj ytterligare set av indata som ska användas för att beräkna koncentrationer baserat på den aktuella standardkurvan.

Dialogrutan **Calculate Additional Concentration** (Beräkna ytterligare koncentrationer) innehåller följande element:

Rullgardinsmenyn Input Data (Indata)	Välj indata för beräkning av ytterligare koncentrationer.
Listan <b>Selected</b> data (Utvalda data)	Listan innehåller namnen på indata för beräkning av ytterligare koncentrationer.
Knappen <b>Add</b> (Lägg till)	Klicka på knappen <b>Add</b> (Lägg till) för att lägga till de aktuellt valda indata i rullgardinsmenyn <b>Input data</b> (Indata) till listan <b>Selected data</b> (Utvalda data).
Knappen <b>Remove</b> (Ta bort)	Klicka på knappen <b>Remove</b> (Ta bort) för att ta bort de aktuellt utvalda data från listan <b>Selected data</b> (Utvalda data).

## Fliken Analysis Type (Analystyp)

Använd den här fliken för att välja analystyp. Se kapitel 11.4 Typer av analys av standardkurvor för en detaljerad beskrivning av analystyperna.

Alternativknappen Analysis type	Välj vilken algoritm för interpolering som ska användas vid beräkning av standardkurvan:
(Analystyp)	Alla analystyper listas nedan och kan väljas:
	1. Point to point (Punkt-till-punkt)
	2. Linear Regression (Linjär regression)
	3. Non-linear Regression (Icke-lineär regression)
	4. Cubic spline (Kubisk spline)
	5. Akima (Akima spline)

	<ol> <li>Polynomial (Polynom) (inklusive urvalsfältet för polynomordning och inklusive viktalternativ)</li> </ol>
	<ol> <li>Four Parameters (Fyra parametrar) (inklusive knappen mer för fastställande av minimum och maximum)</li> </ol>
	<ol> <li>Four Parameters Marquardt (Fyra parametrar Marquardt) (inklusive knappen mer för att definiera viktningsalternativen).</li> </ol>
	<ol> <li>Five parameters (Fem parametrar) (inklusive knappen mer för att definiera viktningsalternativen).</li> </ol>
	10.LogitLog (inklusive knappen mer för fastställande av minimum och maximum).
Rullgardinsmenyn Data scaling (Dataskalning)	Analystypen tillämpas på skalade värden. Följande skalningslägen kan väljas:
	<ul> <li>Lin(x)Lin(y): x- och y-axlarna skalas linjärt</li> </ul>
	<ul> <li>Lin(x)Log(y): x-axeln skalas linjärt, y-axeln skalas logaritmiskt</li> </ul>
	<ul> <li>Lin(x)Log(y): x-axeln skalas logaritmiskt, y-axeln skalas linjärt</li> </ul>
	<ul> <li>Log(x)Log(y): x- och y-axlarna skalas logaritmiskt</li> </ul>

i

# Anmärkning

Observera att om man väljer logaritmisk skalning för x-axeln är det inte möjligt att ha ett koncentrationsvärde på 0, eftersom det matematiskt sett inte är möjligt att beräkna en logaritm av 0. Det är dock möjligt att använda ett värde på till exempel 0.0000001.

Knappen <b>More</b> (Mer)	Klicka på knappen <b>More (Mer)</b> för att visa följande alternativ:
	<ul> <li>Numeriska urvalsfältet Standard Curve (Standardkurva)</li> <li>Om det finns fler än en försöksgrupp för LogitLog och Four Parameters (Fyra parametrar), kan Min och Max väljas individuellt för varje kurva.</li> </ul>
	<ul> <li>Weighting (Viktning)</li> <li>Välj Use weights (Använd vikter) för att använda en av följande viktningsmetoder:</li> </ul>
	<ul> <li>Automatisk med varians</li> <li>Automatisk med relativ vikt</li> <li>Manuell</li> </ul>
	Se kapitel 11.4.14 Viktning för fyra/fem parametrar – Marquardt/Polynomial Fit
	<ul> <li>Dialogrutan Viktning – välj knappen More (Mer) om "Automatic using variance" (Automatisk med varians) är vald=&gt; Dialogrutan Felhantering visas – Definiera de viktningsfaktorer som ska användas om ett av de två fellägena (Alla replikat är lika eller endast ett replikat är kvar) inträffar, vilket ger en varians på 0 och därför skulle leda till ett beräkningsfel.</li> </ul>
	<ul> <li>Textfälten Min/Max Min/Max gör det möjligt för användaren att definiera minsta eller maximala gränsen för standardkurvan för LogltLog och Four Parameters (Fyra parametrar).</li> </ul>



Kryssrutan Include (0,0) (Inkludera (0,0))	Om alternativet <b>Include (0,0) (Inkludera (0,0))</b> är valt läggs (0,0) till som standardpunkt. Detta är endast tillgängligt om <b>Linear (Linjär)</b> har valts som axeldelning för båda axlarna.
Kryssrutan Extrapolation (Extrapolering)	Extrapolering tillämpar koncentrationsberäkningarna på de baspunkter som ligger utanför det tillåtna intervallet samt på de som ligger inom intervallet, till exempel: en extrapoleringsfaktor på 3 tvingar en koncentrationsberäkning av värden mellan min-2*(max- min) och max+2*(max-min), där min och max motsvarar det minsta och maximala koncentrationsvärdet i standardkurvan.
Numeriska fältet Extrapolation factor (Extrapolerings- faktor)	Datafältet är endast aktivt om kryssrutan <b>Extrapolation</b> factor (Extrapoleringsfaktor) har markerats. Den definierar de nya gränserna för koncentrationsberäkningen.

## Fliken Intercepts (Skärningspunkter)

Använd den här fliken för att beräkna koncentrationer för valda Y-värden baserat på standardkurvan.

Spinnkontrollen Exp. group (Försöksgrupp)	Om det finns flera försöksgrupper kan man välja en lämplig grupp.
Rullgardinsmenyn Input Data (Indata)	Alla data som är tillgängliga för beräkning visas i denna lista.
Listfältet Intercept name and formula (Namn på skärningspunkt och formel)	Ett namn på skärningspunkt måste först skapas. Därefter måste formeln för skärningspunkten definieras. Ett numeriskt värde eller en formel kan anges.
Operatorerna Formula input (Inmatning av formel)	<ul> <li>Urvalsrutorna gör det enkelt att mata in formler genom att tillhandahålla en mängd olika funktioner, operatorer och variabler.</li> <li>Variables (Variabler) Alla variabler som accepteras av programmet är tillgängliga i listrutan.</li> <li>Knappen Klicka på den här knappen för att öppna dialogrutan Define Constants (Definiera konstanter).</li> </ul>
	<ul> <li>Operators (Operatorer) Alla operatorer som accepteras av programmet visas här.</li> <li>Functions (Funktioner) Alla funktioner som accepteras av programmet visas här.</li> </ul>

#### Exempel

#### IC50 för standardkurvan:

Formel: (ST1\_1+ST1\_8)/2

värdet av den minsta standarden plus värdet av den högsta standarden, dividerat med 2.



# Fliken Axis (Axel)

Använd den här fliken för att definiera hur axlarna ska se ut.

Grupprutan X-axis (X-axel) Grupprutan Y-axis (Y-axel)	<ul> <li>Följande element finns tillgängliga i varje grupp:</li> <li>Textfältet Label (Etikett): Ange en axeletikett</li> <li>Knappen Color (Färg): Välj en färg för axelns alla element.</li> </ul>
	<ul> <li>Kryssrutan Log-scaling (Logaritmisk skalning) Välj den här kryssrutan för att visa diagrammet axlar logaritmiskt (påverkar inte beräkningen).</li> </ul>
	<ul> <li>Alternativknappen Auto select range (Välj intervall automatiskt): Programvaran fastställer automatiskt minsta och maximala värde.</li> </ul>
	<ul> <li>Alternativknappen Range (Intervall): Numeriska fälten Min och Max är aktiverade. Det minsta och maximala värdet för axeln kan anges.</li> </ul>
	<ul> <li>Kryssrutan Grid (Rutnät): Om detta väljs visas rutnätet på axeln. Knappen Color (Färg) och rullgardinsmenyn Line style (Linjeformat) kan användas för att anpassa rutnätet.</li> </ul>

# Fliken Graph (Diagram)

Använd den här fliken för att definiera diagrammets utseende.

Grupprutan Title (Titel)	<ul> <li>Följande element är tillgängliga:</li> <li>Textfältet Label (Etikett): Ange en etikett för diagrammet.</li> <li>Knappen Color (Färg): Välj en färg för diagrammets etikett.</li> </ul>
Grupprutan <b>Curves (Kurvor)</b>	<ul> <li>Följande element är tillgängliga:</li> <li>Spinnkontrollen Exp. group (Försöksgrupp) Välj den försöksgrupp som ska anpassas (endast tillgänglig om fler än en försöksgrupp har definierats).</li> <li>Knappen Color (Färg): Välj en färg för kurvan för den valda försöksgruppen.</li> <li>Rullgardinsmenyn Symbol: Välj en symbol för baspunkterna i den valda försöksgruppen.</li> <li>Textfältet Label (Etikett): Ange en etikett för kurvan för den valda försöksgruppen.</li> <li>Kryssrutan Hide curve (Dölj kurva) Om den är vald visas inte kurvan.</li> <li>Spinnkontrollen Line Width (Linjebredd): Välj linjebredd för den valda försöksgruppen.</li> </ul>
Grupprutan Font (Teckensnitt)	Välj Small (Liten), Medium (Mellan) eller Large (Stor) teckenstorlek.

Grupprutan <b>Display (Visa)</b>	Följande element kan väljas för att visas: Legend (Förklaring): Kurvfärg, stil för baspunkt och etikett visas tillsammans med ytterligare information (parametrar för analystyp, korrelationskoefficient och så vidare).
	Base points (Baspunkter)
	Intercepts (Skärningspunkter): Om skärningspunkter

anges kan de visas och märkas i diagrammet.

**Error bars (Felstaplar)**: Om standarder definieras som replikat visas en stapel som visar intervallet plus/minus standardavvikelse för varje baspunkt.

# 4.3.9 Concentration transformation (Koncentrationstransformationer): Lägg till nya koncentrationstransformationer

# I dialogrutan Concentration transformations

**(Koncentrationstransformationer)** kan en koncentrationsformel definieras. Formeln används för att transformera indata för koncentration individuellt för varje brunn.

#### I dialogrutan Concentration transformations

(Koncentrationstransformationer) kan ytterligare beräkningar utföras på indata för koncentration.

Fönsterelementen är jämförbara med indata för **transformation** Se kapitel 4.3.5 Transformed data (Transformerad data): Lägg till ny transformation för mer information.

# 4.3.10 Utvärdera data: Definition av cutoff

Använd det här alternativet för att kategorisera antingen rådata eller beräknade data vid tröskelgränser. Tröskelvärden kan definieras som fasta numeriska värden eller formler.

Använd fönstret **Cutoff Definition (Definition av cutoff)** för att definiera cutoffintervall och för att tilldela färger och namn till cutoff-resultaten.

Fönstret Cutoff Definition (Definition av cutoff) innehåller följande element:

Rullgardinsmenyn Input data (Indata)	Välj den Input Data (Indata) som ska användas för utvärderingen.
Urvalslistan <b>Exp. group</b> (Försöksgrupp)	Om plattan innehåller flera tester ska den relevanta försöksgruppen där cutoff ska gälla, väljas ut. Om plattan endast innehåller ett test är endast en grupp tillgänglig och fältet är inte synligt.

I listan **Cutoff** kan upp till tio cutoff-intervall definieras. Gränserna för varje intervall kan anges.

Det högsta gränsvärdet måste anges överst i listan. Den nästa lägsta placeras därefter och så vidare. En lågt-högt-pil visar nivån. Gränsvärdena hör till de övre gränserna (större och lika villkor).

Knapparna <b>Colors</b> (Färger)	En färg kan tilldelas de individuella resultatnivåerna. Vid visning av resultaten kommer dessa färger att ingå och representera de kvalitativa resultaten.
Textfälten	Resultatindikatorerna för varje värdeintervall måste
<b>Labels</b> (Etiketter)	namnges, t.ex.: positiv (pos), negativ (neg), gränsfall(?),

Textfälten Limits (Gränser)	Gränserna kan anges som ett konstant värde eller som en formel. Formler kan skrivas in direkt i ett fält eller via fälten <b>Formula input (Inmatning av formel)</b> . Maximalt nio gränser får definieras.
Grupprutan <b>Formula input</b> (Inmatning av formel)	<ul> <li>Formler som ska skrivas in i fältet Limits (Gränser) kan sammansättas med hjälp av variablerna, operatorerna och funktionerna nedan. Se kapitel 11.3 Hur man skriver en formel för ytterligare detaljer.</li> <li>Rullgardinsmenyn Variables (Variabler) Alla tillgängliga variabler visas.</li> <li>Knappen Klicka på den här knappen för att öppna dialogrutan Define Constants (Definiera konstanter).</li> <li>Rullgardinsmenyn Operators (Operatorer) Alla tillgängliga operatorer visas.</li> <li>Rullgardinsmenyn Functions (Funktioner) Följande funktioner kan väljas.</li> </ul>
Kryssrutan <b>Competitive Test</b> (Kompetitivt test)	I kompetitiva tester tilldelas ett positivt resultat låga värden och ett negativt resultat höga värden. I utvärderingen kommer gränserna att löpa uppifrån och ner och motsvarande resultatsymbol kommer att tilldelas när ett värde är lika med eller mindre än gränsvärdet. Med tanke på detta kommer <b>lågt-högt-pilen</b> att vändas om.
Knappen Cutoff results selection (Urvalet cutoff- resultat)	Klicka på knappen <b>Cutoff results selection</b> (Urvalet cutoff-resultat) för att öppna dialogrutan <b>Cutoff results selection</b> (Urvalet cutoff-resultat).

# Cutoff results selection (Urvalet cutoff-resultat)

Använd det här fönstret för att välja om ett kvalitativt resultat för en viss identifieringstyp ska visas eller inte.

Cutoff results selection... (Urvalet cutoff-resultat) innehåller följande delar:

Trädstrukturen <b>Cutoff display</b> selection (Urvalet cutoff- display)	Alla definierade identifieringsnamn visas i en trädstruktur, grupperade efter identifieringstyper. Varje identifierare är associerad med en kryssruta.
	Om du markera eller avmarkerar kryssrutan för ett identifieringsnamn anger du om de kvalitativa resultaten av de aktuella brunnarna kommer att visas eller inte. Efter utvärderingen visas endast cutoff-resultat för de valda identifieringstyperna.

# 4.3.11 Utvärdera data: QC-validering

Valideringar används för att kontrollera giltigheten hos ett test. Om de definierade kriterierna inte uppfylls visas ett felmeddelande efter mätningen och inga cutoffresultat visas. (Förutom att användaren har rätt att **fortsätta utvärderingen vid felmeddelande**) – se kapitel 9.6 Användarrättigheter.

När du har valt rätt indata ska de önskade formlerna anges i listfältet **Validation Conditions (Valideringsvillkor)**. En formelredigerare hjälper användaren att skapa motsvarande matematiska formler.

#### Exempel:

Om valideringskriterierna definieras på följande sätt:

#### NC1>0 och NC1<0.1

kontrollerar utvärderingen om den negativa kontrollen kan hittas inom det givna intervallet. Om inte visas ett felmeddelande.

Dialogrutan **Define QC Validations (Definiera QC-valideringar)** innehåller följande element:

Rullgardinsmenyn Input Data (Indata)	Välj de relevanta data som valideringen ska tillämpas. Till exempel: Rådata, osv.
Urvalsfältet <b>Exp. group</b> (Försöksgrupp)	Om plattan innehåller flera tester ska den relevanta försöksgruppen där valideringen ska gälla, väljas ut. Om plattan endast innehåller ett test är endast en grupp tillgänglig.
Urvalsfältet Validation group (Valideringsgrupp)	Använd <b>Validation groups</b> ((Valideringsgrupp) för att definiera valideringskriterier för samma försöksgrupp med olika indata.
Listfältet Validation Conditions (Valideringsvillkor)	Denna endimensionella lista kommer att fyllas med de formler och logiska ekvationer som definierar <b>Validation</b> <b>Conditions</b> (Valideringsvillkor). Dessa logiska ekvationer ger ett logiskt resultat. Programmet undersöker indata med hjälp av denna ekvation, och om valideringskriterierna är uppfyllda ger programmet resultatet TRUE. Om kriterierna inte uppfylldes visas resultatet FALSE och ett felmeddelande.
Grupprutan <b>Formula input</b> (Inmatning av formel)	<ul> <li>Delar av formlerna som ska anges i fältet Validation Conditions (Valideringsvillkor) kan väljas från följande rullgardinsmenyer:</li> <li>Rullgardinsmenyn Variables (Variabler) Alla tillgängliga variabler visas.</li> <li>Knappen Klicka på den här knappen för att öppna dialogrutan Define Constants (Definiera konstanter).</li> <li>Rullgardinsmenyn Operators (Operatorer) Alla tillgängliga operatorer visas.</li> <li>Rullgardinsmenyn Functions (Funktioner) Med hjälp av den angivna paletten, kan följande funktioner väljas:</li> <li>Se kapitel 11.3 Hur man skriver en formel för ytterligare detaljer.</li> </ul>



# Plate to Plate QC (QC platta-till-platta)

Använd det här alternativet för att definiera en långsiktig QC-validering (se kapitel 7.4.2 Verktygsfältet: File (Fil)/QC platta-till-platta):

Rullgardinsmenyn Input data (Indata)	Välj de relevanta data som valideringen ska tillämpas från rullgardinsmenyn.
Kryssrutor <b>Kontroll</b>	Välj typ av kontroll från rullgardinsmenyn och ange sedan det förväntade medelvärdet och standardavvikelsen.
	En rullningslist visas bredvid dialogkontrollerna om du vill definiera fler än fyra QC-kontroller.
Alternativen Workspaces to be evaluated (Arbetsytor som ska utvärderas)	Välj lämpligt filter för att definiera vilka arbetsytor som ska utvärderas.

# 4.3.12 Datahantering: Dataexport

Data som ska exporteras till en ASCII- eller Excel-fil kan väljas i den här dialogrutan:

Listan <b>Available data</b> (Tillgängliga data)	En lista med alla tillgängliga data visas, beroende på metoddefinitionen.
← och → knappar	Data för export kan väljas enkelt genom att dra och släppa eller genom att klicka på data i fönstret <b>Available</b> <b>data</b> (Tillgängliga data) och sedan på pilen som pekar mot fönstret <b>Selected data</b> (Utvalda data). Data kan avmarkeras genom att utföra proceduren i omvänd ordning
Knapparna <b>Up</b> (Upp) och <b>Down</b> (Ner)	Ordningen på de utvalda data kan ändras genom att välja ett element och trycka på knappen <b>Up</b> (Upp) för att flytta det uppåt eller <b>Down</b> (Ner) för att flytta det nedåt.
Listan <b>Selected data</b> (Utvalda data)	Den här listan visar alla utvalda data som har överförts från fältlistan <b>Available data</b> (Tillgängliga data).
Knappen <b>Export options</b> (Exportera alternativ)	Klicka på den här knappen för att öppna dialogrutan Export options (Exportera alternativ).

Den faktiska exporten av data sker endast med hjälp av automatiserad databehandling (se 4.3.14 Datahantering: Automatiserad datahantering).

# Export Options (Exportera alternativ)

Dialogrutan Export options... (Exportera alternativ) innehåller följande element

Grupprutan Direction (Riktning)	Användaren kan bestämma om plattdata ska extraheras horisontellt rad för rad eller vertikalt, kolumn för kolumn, och skrivas i denna ordning till filen.
	<ul> <li>Alternativknappen Horizontal (Horisontell) Data samlas i rader.</li> </ul>
	<ul> <li>Alternativknappen Vertical (Vertikal) Data samlas i kolumner.</li> </ul>
Grupprutan Result (Resultat)	Välj om du vill exportera data i form av en matris eller en kontinuerlig lista:
	<ul> <li>Alternativknappen Matrix (nested) (Matris - nästlad) Data från alla utvalda dataset ordnas i en enstaka matris. Denna matris innehåller de första kolumnerna för alla dataset följt av de andra kolumnerna för alla dataset.</li> </ul>
	<ul> <li>Alternativknappen Matrix (separated) (Matris - separerad)</li> <li>Data för varje utvald dataset ordnas i en separat matris.</li> </ul>
	<ul> <li>Alternativknappen Matrix (XFluor style) (Matris - Xfluor-stil)</li> <li>Data för varje utvald dataset ordnas i en separat matris tillsammans med kolumn- och brunnsbeskrivning på liknande sätt som i XFluor.</li> </ul>
	<ul> <li>Alternativknappen Table (well data in rows) (Tabell - brunnsdata i rader)</li> <li>Alla data för en brunn ordnas i horisontell riktning, och börjar med data i brunnen som är placerad i A1.</li> </ul>
	<ul> <li>Alternativknappen Table (well data in columns) (Tabell - brunnsdata i kolumner) Alla data för en brunn ordnas i vertikal riktning, och börjar med data i brunnen som är placerad i A1.</li> </ul>
	Om kryssrutan <b>Add kinetic timestamps (Lägg till kinetiska tidsstämplar)</b> är markerad läggs tidsstämplarna för mätningarna till.
	Om kryssrutan <b>Add Temperatures (Lägg till</b> temperaturer) är markerad läggs temperaturen för mätningarna till.
	Om kryssrutan <b>Insert Data names (Infoga datanamn)</b> är markerad visas en lista med namnen på exporterade data ovanpå ASCII-filen respektive Excel-kalkylbladet.


Grupprutan Add data (Lägg till data)	<ul> <li>Ger valbara dataalternativ.</li> <li>Den valda informationen bifogas i slutet av data och exporteras sedan tillsammans med utvalda data.</li> <li>Välj respektive kryssrutor:</li> <li>Kryssrutan Date/time of measurement (Datum/tid för mätning)</li> <li>Kryssrutan Method filename (Filnamn på metod)</li> <li>Kryssrutan Method pathname (Sökvägsnamn på metod)</li> <li>Kryssrutan Workspace filename (Filnamn på arbetsyta)</li> <li>Kryssrutan Workspace pathname (Sökvägsnamn på arbetsyta)</li> <li>Kryssrutan Filter wavelength value(s) (Filtrera våglängdsvärden)</li> <li>Kryssrutan Current user name (Nuvarande användarnamn)</li> <li>Kryssrutan Multiplate plate information (Information om flera plattor)</li> <li>Kryssrutan Workspace audit trail (Arbetsyta audit trail) (endast magellan Tracker)</li> <li>Kryssrutan Workspace signatures (Underskrifter i arbetsytor) (endast magellan Tracker)</li> </ul>
Knappen Set as default (Ställ in som standard)	Inställningarna kan sättas som standardinställningar för framtida användning.
Restore default button (Knappen Återställ till standard)	Inställningarna återställas till de tidigare definierade standardvärdena.
Exempel	
Rådata, kinetikcykel 1	, tidsstämpel 0 s
11 12 13	
21 22 23	
Rådata, kinetikcykel 2	2, tidsstämpel 33 s
81 82 83	
91 92 93	

Matris (nästlad), horisontell: 11 81 12 82 13 83 21 91 22 92 23 93	Matris (separerad), horisontell, med tidsstämplar: 0 s 11 12 13 21 22 23 33 s 81 82 83 91 92 93	Matris (XFluor-stil): <> 1 2 3 A 11 12 13 B 21 22 23 <> 1 2 3 A 81 82 83 B 81 82 83
Tabell (brunnsdata i rader), horisontell, med tidsstämplar: 0 s 33 s 11 81 12 82 13 83 21 91 22 92 23 93	Tabell (brunnsdata i kolumner), horisontell: 11 12 13 21 22 23 81 82 83 91 92 93	Tabell (brunnsdata kolumner), vertikal: 11 21 12 22 13 23 81 91 82 92 83 93

# Export to ASCII File (Exportera till ASCII-fil)

Dialogrutan Export to ASCII File (Exportera till ASCII-fil) innehåller följande element:

Grupprutan <b>Decimal character</b> (Decimaltecken)	<ul> <li>Decimaltecknet kan vara en punkt eller ett kommatecken:</li> <li>Alternativknappen Point (x.xx) (Punkt)</li> <li>Alternativknappen Comma (x,xx) (Komma)</li> </ul>
Grupprutan <b>Delimiter</b> (Avgränsare)	Det här alternativet ställer in en avgränsare för att separera de individuella värdena inom data på raderna i en ASCII-fil:
	<ul> <li>Alternativknappen <b>Tabulator</b> (Tabulator, tabb) Standardalternativet använder tabb som avgränsare.</li> </ul>
	<ul> <li>Alternativknappen Other symbol (Annan symbol) Om det väljs kan ett användardefinierat avgränsningstecken anges.</li> </ul>
Grupprutan <b>Path</b> (Sökväg)	<ul> <li>Det här alternativet anger sökvägen där ASCII-filen kommer att lagras:</li> <li>Alternativknappen Export default (Exportera standard)</li> <li>Alternativknappen Use path (Använd sökväg): Välj en sökväg där ASCII-filen ska lagras.</li> </ul>
Grupprutan <b>Encoding</b> (Kodning)	<ul> <li>Välj exportkod:</li> <li>ANSI: För ANSI-kompatibla tecken.</li> <li>Unicode: För tecken som inte kan exporteras med ANSI-kod (t.ex. kinesiska, kyrilliska).</li> </ul>



Knappen Set as default (Ställ in som standard)	Inställningarna kan sättas som standardinställningar för framtida användning.
Knappen <b>Restore default</b> (Återställ till standard)	Inställningarna kan återställas till de tidigare definierade standardvärdena.



# FÖRSIKTIGHET

EN LÄMPLIG AVGRÄNSARE FÖR EXPORT MÅSTE VÄLJAS. EXPORTAVGRÄNSAREN OCH DECIMALTECKNET FÅR INTE VARA DETSAMMA. TAB ÄR STANDARD SOM AVGRÄNSARE, MEN EN ANNAN SYMBOL KAN VÄLJAS.

# Export to Excel (Exportera till Excel)

Dialogrutan Export to Excel (Exportera till Excel) innehåller följande element:

Grupprutan <b>Target</b> (Mål)	Alternativknapparna definierar var överförda data ska placeras i Excel. Detta alternativ gäller även för automatisk Excel-export.
	<ul> <li>Alternativknappen New workbook (Ny arbetsbok) Överförda data lagras i det första kalkylbladet i en ny arbetsbok. En ny fil skapas alltid, oavsett om Excel startas eller inte.</li> </ul>
	<ul> <li>Alternativknappen New worksheet (Nytt kalkylblad) De överförda data placeras i ett nytt kalkylblad i en öppen, aktiv Excel-arbetsbok. Om Excel inte är öppet när du väljer det här alternativet startas det och en ny arbetsbok skapas med önskade data i det första kalkylbladet.</li> </ul>
	<ul> <li>Alternativknappen Insert into worksheet at cell (Infoga i kalkylbladet i cell) och textfält för cellkoordinater</li> <li>Det första värdet av överförda data placeras i den angivna cellen (standard är cell A1) i ett öppet, aktivt Excel-kalkylblad.</li> </ul>
	Om Excel inte är öppet när du väljer det här alternativet kommer det att startas och en ny arbetsbok skapas. Önskade data infogas i det första kalkylbladet i den definierade cellen.
	<ul> <li>Alternativknappen Append to current worksheet (Bifoga till aktuellt kalkylblad) Med det här alternativet bifogas data till det aktuella kalkylbladet. Om Excel inte är öppet när du väljer det här alternativet kommer det att startas och en ny arbetsbok skapas. Önskade data infogas i det första kalkylbladet.</li> </ul>
	<ul> <li>Använd Insert into Template (Infoga i mall) för att exportera data till en fördefinierad Excel-mall. Sökväg och namn på Excel-mallen måste anges genom att klicka på knappen Motsvarande cellposition måste anges.</li> </ul>

Knappen Set as default (Ställ in som standard)	Inställningarna kan sättas som standardinställningar för framtida användning.
Knappen <b>Restore default</b> (Återställ till standard)	Inställningarna kan återställas till de tidigare definierade standardvärdena.

### 4.3.13 Datahantering: Utskriven rapport

Dialogrutan Printed report (Utskriven rapport) innehåller formateringsfunktioner för pappersutskrifter för att anpassa innehållet och utseendet på utskriften.

Inställningarna för rapporterna är grupperade i fyra flikar som representerar följande formateringsområden:

- Fliken Data Selection (Dataurval)
- Fliken Page Setup (Sidinställning)
- Fliken Header (Sidhuvud)
- Fliken Footer (Sidfot)

#### Fliken Data Selection (Dataurval)

På fliken Data Selection (Dataurval) finns en lista över data som kan skrivas ut.

Skriv ut som gruppruta	Välj om du vill skriva ut data i form av en matris eller en lista:
	<ul> <li>List (Lista) Det individuella listformatet (horisontellt, vertikalt, sorterat och alternerande) måste konfigureras i dialogrutan Properties (Egenskaper) (se nedan).</li> </ul>
	<ul> <li>Matrix (Matris) Med matrisvalet skrivs data ut i en tvådimensionell lista (matris).</li> </ul>
Lista över datakällor och destinationer	<ul> <li>Två listor används för att definiera de utskrivna data:</li> <li>Listan Available data (Tillgängliga data) Lista över alla tillgängliga data som kan skrivas ut. Sidbrytningar och skiljelinjer kan läggas till för att uppnå önskad layout. Element som ska skrivas ut kan väljas genom att dubbelklicka; de finns i listan över valda data.</li> </ul>
	<ul> <li>Listan Selected data (Utvalda data)         Lisa över all utvalda data som kommer att skrivas ut.         Informativa anteckningar hjälper också till att förklara         hur data kommer att skrivas ut, t.ex. om det kommer             att visas i form av en matris, en lista, som             datainsamling eller som ett diagram. Poster som visas             under en rubrik kommer att inkorporeras i             motsvarande list- eller matrisformat, vilket gör det             möjligt att jämföra olika sets av värden, t.ex.             mätvärden och resultat.         </li> </ul>
Knappen <b>Append</b> (Bifoga)	Klicka på den här knappen för att överföra data som markerats i fönstret <b>Available data (Tillgängliga data)</b> till fönstret <b>Selected data (Utvalda data)</b> . Du kan också använda <b>drag and drop (Dra och släpp)</b> .



# 4. Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)

Knappen Insert (Infoga)	Använd det här alternativet om du vill visa två set av likadant formaterade data inom samma område. Först måste data från listan <b>Available data (Tillgängliga data)</b> väljas, sedan data från listan <b>Selected data (Utvalda data)</b> . När du klickar på knappen <b>Insert (Infoga)</b> kommer det förstnämnda elementet att inkluderas i det sistnämnda och det kommer då att uppta samma område i utskriften som det överordnade elementet. Om t.ex. standardavvikelsen ingår i elementets mätvärde, visas båda i den motsvarande cellen av en matris. Mätvärdena kommer i detta fall att fungera som det överordnade elementet och kommer därför att visas först.
Knappen <b>Remove</b> (Ta bort)	Valda element i listan <b>Selected data (Utvalda data)</b> kan raderas.
Knapparna <b>Up</b> (Upp) och <b>Down</b> (Ner)	Använd dessa knappar för att omorganisera sekvensen av listor, diagram, matriser på utskriften eller för att omorganisera elementet inom en matris eller en lista.
Knappen <b>Properties</b> (Egenskaper)	Aktiveras när en lista väljs. Du kan välja stil och innehåll i listan.
Gruppruta <b>Style</b> (Stil)	Alternativknappen <b>Vertical list (Vertikal lista)</b> Skriver ut data i form av en vertikal lista (ordning A1, B1 och så vidare).
	Alternativknappen Vertical list alternating (Vertikal lista alternerar) Ordningen av data alternerar för varje rad eller kolumn (ordning A1, B1, H1, H2, G2, A2)
	Alternativknappen <b>Horizontal list (Horisontell lista)</b> Skriver ut data i form av en horisontell lista (ordning A1, A2 och så vidare).
	Alternativknappen Horizontal list alternating (Horisontell lista alternerar) Ordningen av data är alternerande för varje rad eller kolumn (ordning A1, A2,A12, B12, B11,B1).
	Alternativknappen <b>Sorted list (Sorterad lista)</b> Det här alternativet sorterar poster i alla kolumner enligt posterna i den första kolumnen. De alfanumeriska posterna i den första kolumnen är uppdelade i två delar, varav den första innehåller endast bokstäver och den andra siffror. Poster sorteras först i alfabetisk ordning. Poster med samma namn sorteras sedan numeriskt. Exempel: Du kan använda det här alternativet med prov- ID:n eller med layouten, begränsat till en försöksgrupp. En sortering av numeriska värden fungerar endast om alla poster är heltal.



# Knappen Sample oriented list (Orienterad lista för prov)

Gör det möjligt att skriva ut individuella listor per prov. Separationen för de enstaka tabellerna görs antingen med hjälp av prov-ID:n (om tillgängligt) eller med hjälp av identifieringsnamnen på plattlayouten. Prov-ID:n behövs om layouten innehåller olika försöksgrupper som representerar olika parametrar för samma prover. Det är möjligt att kombinera listformatet **Sample oriented list (Orienterad lista för prov)** med andra **parametrar** som identifierare, försöksgrupper och cutoff-resultat.

yle vamole				Limitations	OK
Layo Raw Blaz Cub 1 A1 A2 Layo Raw	out data bit results 2 BL1_1 0.1 BL1_2 0.1	3 4 01~ -0.000 02~ 0.000	<ul> <li>Vertical list</li> <li>Vertical list alternating</li> <li>Horizontal list</li> <li>Horizontal list alternating</li> <li>Sorted list</li> <li>Sample oriented list(s)</li> </ul>	Identifiers         SM         BL         ST         PC         Exp. Group(s)         Image: All experimental groups	Cance Help
Cub	aff results			🔾 0 nly 👔 🗸 🗸	
1	2	3			
A3 A4	SM1_2_1 SM1_2_3	0.103~ 0.104~		Cutoff results	
B3 B4	SM1_2_2 SM1_2_4	0.203 0.204		O All Cutoff results	
			4		

Grupprutan	Välj mellan följande alternativ:
Limitations (Begränsningar)	Identifiers (identifierare): välj eller avmarkera vilka identifierare som ska visas i listan.
	<b>Exp. Group (Försöksgrupp)</b> : välj mellan alla försöksgrupper eller ange numret på den försöksgrupp som ska visas i listan.
	Cutoff results (Cutoff-resultat): välj mellan alla cutoff- resultat eller välj från listan vilka resultat som ska visas i

listan.



# Fliken Page Setup (Sidinställning)

På fliken **Page Setup (Sidinställning)** definieras allmänna inställningar:

Grupprutorna Display Header/Footer (Visa sidhuvud/sidfot)	<ul> <li>Kryssrutan Print Header (Skriv ut sidhuvud) Om det väljs, skrivs sidhuvudet ut.</li> <li>Kryssrutan Print Footer (Skriv ut sidfot) Om det är valt, skrivs sidfoten ut.</li> <li>Alternativknappen Header (Sidhuvud) <ul> <li>Every page (Varje sida) (skriver ut sidhuvudet på varje sida)</li> <li>First page only (Endast första sidan) (skriver ut sidhuvudet endast på den första sidan)</li> <li>Except first page (Förutom första sidan) (skriver ut sidhuvudet på alla sidor utom den första)</li> </ul> </li> <li>Alternativknappen Footer (Sidfot) <ul> <li>Every page (Varje sida) (skriver ut sidfoten på varje sida)</li> <li>First page only (Endast första sidan) (skriver ut sidfoten på alla sidor utom den första)</li> </ul> </li> </ul>
Gruppen Print/Page (Skriv ut/sida)	<ul> <li>Alternativknappen Color (Färg) Utskriften är i färger, om skrivaren har stöd för det. Brunnsspecifika data skrivs ut i layoutens färg (till exempel: rådata från provet skrivs ut i blått). Cutoff- resultaten skrivs ut i intervallets färg (till exempel: "pos"-brunnar skrivs ut i rött). Diagrammen skrivs ut så som den visas.</li> <li>Alternativknappen Black and white (Svartvit) Utskriften är svartvit.</li> </ul>
Grupprutan Margins (Marginaler)	<ul> <li>I rullgardinslistan Unit (Enhet) kan du välja cm eller tum. Det finns fyra numeriska fält som definierar marginalerna:</li> <li>Textfältet Left (Vänster): ange värdet för den vänstra marginalen.</li> <li>Textfältet Right (Höger): ange värdet för den högra marginalen.</li> <li>Textfältet Above (Ovanför): ange värdet för den övre marginalen.</li> <li>Textfältet Below (Nedanför): ange värdet för den nedre marginalen.</li> </ul>

# Fliken Header (Sidhuvud)

ata selection	Page setup Header Fr	ooter			
	Option:	Adjustment			Row number:
	🗹 Page numbers	🔿 left	O centered	💿 right	3 🔹
	🗹 Date	<li>left</li>	◯ centered	🔘 right	2 🔹
	🗹 Time	🔘 left	O centered	🔘 right	3
	🗹 File name	Ieft	Centered	🔘 right	1
	Program name	) left	C centered	🔘 right	1
	Comment	Ieft	Centered	🔘 right	1
	Separators				
	Above	Below			

På fliken **Header (Sidhuvud)** definieras innehållet i dokumentsidhuvudet:

Kryssrutan <b>Option (Alternativ)</b>	<ul> <li>Använd Option (Alternativ) för att välja någon av följande kryssrutor som ska skrivas ut i sidhuvudet:</li> <li>Page numbers (Sidnummer)</li> <li>Date (current) (Aktuellt datum)</li> <li>Time (Tid)</li> <li>File Name (Filnamn)</li> <li>Program name (Programnamn)</li> <li>Comment (Kommentar)</li> <li>Textfältet Comment (Kommentar) är aktivt, eftersom kryssrutan Comment (Kommentar) har markerats.</li> </ul>
Alternativknappen Adjustment (Justering)	<ul> <li>Välj ett av följande alternativ:</li> <li>Left (Vänster)</li> <li>Centered (Centrerad)</li> <li>Right (Höger)</li> </ul>
Listan <b>Row number</b> (Radnummer)	Row number (Radnummer) är ett numeriskt fält där raden för varje alternativ kan justeras.
Kryssrutan Separators (Avskiljare)	<ul> <li>I kryssrutorna för avskiljare anges var gränslinjerna ska visas. Dessa kan placeras under, över eller både under och över sidhuvudet.</li> <li>Kryssrutan Above (Ovanför): Avskiljarlinjen visas ovanför rubriken.</li> <li>Kryssrutan Below (Nedanför): Avskiljarlinjen visas under rubriken.</li> </ul>



#### Fliken Footer (Sidfot)

På fliken **Footer (Sidfot)** definieras innehållet i dokumentfoten. Fliken **Footer (Sidfot)** innehåller samma element som fliken **Header (Sidhuvud)**. För en beskrivning av fliken Footer (Sidfot) se fliken sidhuvud.

Fliken Footer (Sidfot) försvinner om **Print Footer (Skriv ut sidfot)** inte är vald på fliken **Page setup (Sidinställning)**.

#### 4.3.14 Datahantering: Automatiserad datahantering

Efter en mätning kan vissa åtgärder startas automatiskt.

Om du väljer dessa åtgärder utförs de kronologiskt i följande ordning:

- Om kryssrutan load sample ID list (Ladda prov-ID-lista) är markerad laddas en prov-ID-lista automatiskt.
- Om kryssrutan save workspace (Spara arbetsyta) är markerad (alltid markerad i magellan Tracker) sparas den skapade arbetsytan automatiskt.
- Om kryssrutan **export to ASCII file (Exportera till ASCII-fil)** är markerad exporteras de data som valts för export automatiskt till en ASCII-fil.
- Om kryssrutan export to ASTM (Exportera till ASTM) är markerad exporteras de data som valts för export automatiskt i ASTM-format efter att metoden körts och kan användas vidare i kompatibla laboratorieinformationssystem (LISLIS - Laboratorieinformationssystem).
- Om kryssrutan **export to Sample Tracking (Export av spårning av prov)** är markerad exporteras de data som valts för export automatiskt till Sample Tracking-systemet (Spårning av prov) när metoden körs. Mer information finns i motsvarande manual för Sample Tracking (Spårning av prov).
- Om kryssrutan **export to Excel (exportera till Excel)** är markerad exporteras de data som valts för export automatiskt till Excel.
- Om kryssrutan **print (Skriv ut)** är markerad skrivs de data som valts för den utskrivna rapporten automatiskt ut på standardskrivaren.
- Om kryssrutan view results after measurement (visa resultat efter mätning) är markerad visas mätdata och beräknade data för granskning.

#### More - Load Sample ID List (Mer - Ladda en prov-ID-lista)

Dialogrutan More (Load Sample ID List) (Mer - Ladda prov-ID-lista) innehåller följande element:

Grupprutan	<ul> <li>Välj filnamnet för den prov-ID-lista som ska laddas:</li> <li>Alternativknappen Same filename as workspace</li></ul>
<b>Filename</b>	(Samma filnamn som för arbetsytan) <li>Alternativknappen Filename (Filnamn) Ange ett</li>
(Filnamn)	filnamn i textfältet.
Grupprutan <b>Path</b> (Sökväg)	<ul> <li>Välj den sökväg där prov-ID-listan lagras:</li> <li>Alternativknappen Sample ID list default path (Standardsökväg för prov-ID-lista)</li> <li>Alternativknappen Use path (Använd sökväg): Välj valfri sökväg.</li> </ul>

#### 4. Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)



Grupprutan Format	Välj formatet för den prov-ID-lista som ska laddas (se kapitel5.2.3 Importera en prov-ID-lista för detaljer):
	<ul> <li>Alternativknappen Autodetect from the list of supported formats (Autodetektera från listan över format som stöds)</li> </ul>
	<ul> <li>Alternativknappen Custom file format (Anpassa filformat): Klicka på knappen File format (Filformat) för att öppna dialogrutan Custom Format (Anpassa format).</li> </ul>



Anmärkning

Om en endast delvis fylld prov-ID-lista laddas automatiskt, raderas alla oanvända prover automatiskt från layouten. Om en slutlig kontroll (t.ex. i brunn H12) har definierats i layouten flyttas denna kontroll automatiskt till brunnen efter den senast använda provbrunnen.



Anmärkning Automatiskt laddade prov-ID-listor bör endast innehålla prov-ID:n och får inte innehålla ID:n för icke-prover (kontroller, standarder osv.).

### More – Workspace Name (Mer – Arbetsytans namn)

När du väljer **Save workspace – More (Spara arbetsyta – Mer)** visas dialogrutan **Workspace Name (Arbetsytans namn)** med följande element:

Grupprutan <b>Format</b>	<ul> <li>En av de föreslagna alternativknapparna måste väljas för att definiera standardnamn för arbetsytan. Namnet kan bestå av datumet, räknaren eller bokstäver. En annan möjlighet är att kombinera datumet med ditt urval av bokstäver. Räknare som används tillsammans med datumet nollställs dagligen, annars ökas räknaren med varje mätning:</li> <li>Datum (DDMMYYYYY) + alternativknappen Counter</li> </ul>
	<ul> <li>Datum (YYYYMMDD) + alternativknappen Counter (Räknare)</li> </ul>
	• Alternativknappen Total Counter (Totalräknare)
	<ul> <li>Alternativknappen Use available barcode (Använd tillgänglig streckkod)</li> </ul>
Textfältet Prefix	Den angivna texten läggs till i början av arbetsytan namn enligt definitionen ovan.
Textfältet <b>Example</b> (Exempel):	Textfältet <b>Example</b> (Exempel) visar ett exempel på det aktuella urvalet.
Knappen <b>Set as default</b> (Ställ in som standard)	Inställningarna kan sättas som standardinställningar för framtida användning.
Knappen <b>Restore default</b> (Återställ till standard)	Inställningarna kan återställas till de tidigare definierade standardvärdena.

# More – Export to LIS (Mer – Exportera till LIS)

Dialogrutan **More (Export to LIS)** (Mer – Exportera till LIS) innehåller följande element:

Grupprutan ASTM delimiter definition (Definition av ASTM- avgränsare)	<ul> <li>Definiera ASTM-avgränsare:</li> <li>Field delimiter (Fältavgränsare): Ange ett tecken för att separera fält (  som standard)</li> <li>Repeat delimiter (Avgränsare för upprepning): Ange ett tecken för upprepning (\ som standard)</li> <li>Component delimiter (Avgränsare för komponent): Ange ett tecken för att separera komponenter (som standard ^).</li> <li>Escape character (Escape-tecken): Ange ett tecken för escape (&amp; som standard).</li> </ul>
Grupprutan <b>Sender ID</b> (Avsändar-ID)	<ul> <li>Definiera inställningarna för Sender ID som ska visas i sidhuvudet på en ASTM-fil:</li> <li>Textfältet Sender ID (Avsändar-ID): Tomt som standard. Om namnet är definierat, t.ex. SUNRISE, visas det i sidhuvudet på ASTM-filen (e.g. H \^&amp;   SUNRISE^magellan V7.2^Tracker).</li> <li>Append app. name (Bifoga programnamn): Om vald exporteras programnamnet, dvs. magellan i sidfoten på ASTM-filen (t.ex. magellan V7.2^Tracker).</li> <li>Append app. version (Bifoga programversion): Om det väljs exporteras versionsnumret av magellan i sidhuvudet på ASTM-filen (t.ex. Standard eller Tracker)</li> </ul>
Grupprutan <b>LIS assay name</b> (LIS-analysnamn)	<ul> <li>Välj det LIS-analysnamn som ska visas i testorder- och resultatposten i ASTM-filen.</li> <li>Use method name (Använd metodnamn): namnet på den metod som används för mätningen.</li> <li>Textfältet Use name (Använd namn): definiera ett nytt namn som ska visas.</li> <li>Exempel: ASTM_Test</li> <li>Testorderpost: O 1 SampleID  ^^^method name  Date and time of measurement (O 1 ProvID   ^^^metodnamn  Datum och tid för mätning)</li> <li>Resultatpost: R 1 ^^ method name^results (R 1 ^^^</li> </ul>
Grupprutan <b>Encoding</b> (Kodning)	<ul> <li>Välj exportkod:</li> <li>ANSI: För ANSI-kompatibla tecken.</li> <li>Unicode: För tecken som inte kan exporteras med ANSI-kod (t.ex. kinesiska, kyrilliska).</li> </ul>
Grupprutan <b>Path</b> (Sökväg)	<ul> <li>Definiera sökvägen för ASTM-export :</li> <li>Export default path (Exportera till standardsökväg): exporten utförs till den mapp som definieras i Miscellaneous (Övrigt).</li> <li>Textfältet Use path (Använd sökväg): definiera en ny exportväg.</li> </ul>



Knappen Set as default (Ställ in som standard)	•	Med det här alternativet kan inställningarna sättas som standardinställningar för framtida användning.
Knappen <b>Restore default</b> (Återställ till standard)	•	Med det här alternativet kan inställningarna återställas till de tidigare definierade standardvärdena.

### ASTM Export – Example File (ASTM-export – Exempelfil)

Varje ASTM-exportfil (.txt) består av följande komponenter:

- Meddelandets rubrikpost: H|\^&|||SUNRISE^magellan V7.2^Tracker
- Patientens informationspost (P): P|1||Sample ID 1 for patient 1 (P|1||Prov-ID 1 för patient 1)
- Testorderpost (O): O|1|SampleID||^^^method name||Date and time of measurement (O|1|ProvID|||^^^metodnamn||Datum och tid för mätning)
- Resultatpost (R): R|1|^^^ method name^results (R|1|^ metodnamn^resultat)
- Meddelandets avslutarpost: L|1|N

#### Exempel:

Inställningar för ASTM-export:

ASTM delimiter definition	Sender ID		OK
Field delimiter:	SUNRISE		Cancel
Repeat delimiter:     \       Component delimiter:     ^       Escape character:     &	Append ap	pp. name pp. version	
LIS assay name		Encoding	
Use method name		<ul> <li>ANSI</li> <li>Unicode</li> </ul>	
Path			
Export default path			
🖲 Use path:			Set as default
C:\Users\Public\Documen	its\Tecan\Magellan\a	isc	Bestore defau



Motsvarande ASTM-fil:

H|\^&|||SUNRISE^magellan V7.2^Tracker

P|1||sample1

O|1|sample1||^^^ASTM\_Test||20180315091504

R|1|^^^ASTM\_Test^Raw data^A1|0.101~|OD|

P|2||sample2

O|1|sample2||^^ASTM\_Test||20180315091504

R|1|^^^ASTM\_Test^Raw data^B1|0.201~|OD|

P|3||sample3

O|1|sample3||^^ASTM\_Test||20180315091504

R|1|^^^ASTM\_Test^Raw data^C1|0.301~|OD|

# More - View Results (Mer - Visa resultat)

Dialogrutan **More – View Results** (Mer – Visa resultat) innehåller följande element:

Grupprutan	Välj de data som ska visas i plattvyn:		
Plate view (Plattvy)	<ul> <li>Rullgardinsmenyn 1st line in well (fixed) (1:a raden i brunnen - fast): Välj ett namn från de tillgängliga datanamnen.</li> </ul>		
	<ul> <li>Rullgardinsmenyn 2nd line in well (fixed) (2:a raden i brunnen - fast): Välj ett namn från de tillgängliga datanamnen.</li> </ul>		
	<ul> <li>Rullgardinsmenyn 3rd line in well (fixed) (3:e raden i brunnen - fast): Välj ett namn från de tillgängliga datanamnen. Till skillnad från de två första raderna tillämpas detta urval endast för första gången. När annan data som ska visas i plattvyn väljs i kontrollfältet visas denna data på 3:e raden. De två första raderna förblir oförändrade.</li> <li>1:a och 2:a raden kan utelämnas genom att välja not</li> </ul>		
	used (Inte använd), vilket gör plattvyn mer lättläst.		
Grupprutan <b>Text box</b> (Textruta)	Välj ett namn från de tillgängliga datanamnen. Detta urval tillämpas endast för första gången. När annan data som ska visas i textrutan väljs i kontrollfältet visas denna data på 3:e raden.		

# 4.3.15 Miscellaneous (Övrigt): Användarprompt

# l Kontrollfältet, expandera Miscellaneous (Övrigt) och välj User prompts (Användarprompt).

l varje **Prompt**-fält i listan kan du ange ett nyckelord. Nyckelordet visas sedan före mätningen. Text kan skrivas in, t.ex. LOT-nummer eller analytikerns fullständiga namn.

Genom att välja **Req. (required) (Krav)** kan en mätning endast påbörjas efter mottagande av en inmatning (en textinformation måste läggas till i fältet **Prompt answer (Promptsvar)**).

Prompter kommer också att visas i listan över tillgängliga data för utskrift.

# 4.3.16 Miscellaneous (Övrigt): Nummerformat

I Kontrollfältet, expandera Miscellaneous (Övrigt) och välj Number format (Nummerformat).

Dialogrutan **Number format (Nummerformat)** används för att ställa in det numeriska formatet för visning av värden. Dessutom kan det vetenskapliga talformatet för värden utanför intervallet väljas.

Standardinställningen är 5 signifikanta siffror. Om du väljer **User defined** (Användardefinierad) kan du ställa in ett sifferformat med 1–10 heltalsställen och 0–7 decimaler.

Om det användardefinierade nummerformatet har valts måste de två numeriska datafälten ställas in.

Dessutom kan den vetenskapliga (med exponent) eller icke-vetenskapliga beskrivningen användas. Ett exempel på formatet för det aktuella valet visas.

Klicka på Set as default (Ställ in som standard) eller Restore default (Återställ till standard) för att spara den inmatade definitionen.

## 4.3.17 Miscellaneous (Övrigt): Metodanteckningar

l Kontrollfältet, expandera Miscellaneous (Övrigt) och välj Method notes (Metodanteckningar).

I fältet **Method notes (Metodanteckningar)** kan användaren lägga till några kommentarer till filnamnet som visas på metodutskriften.

Enkel formatering som fet, kursiv, understruken och färgad text är möjlig. Formatering kan utelämnas vid utskrift.

# 4.4 Spara metoden

Klicka på Finish (Slutför) och spara metoden via fönstret Save as (Spara som), som innehåller följande element:

Fältet <b>Filename</b> (Filnamn)	Ange eller ändra metodens filnamn.
Fältet File remarks (Filkommentarer)	Lägg till några kommentarer till filnamnet, som kommer att visas till exempel i dialogrutorna <b>Open file</b> (Öppna fil).
Fältet <b>Audit trail comment</b> (Audit trail kommentar)	Lägg till en kommentar, som kommer att sparas i audit trail (endast tillgängliga för <b>magellan</b> Tracker).
Knappen Organize favorites (Organisera favoriter)	Klicka på den här knappen för att öppna dialogrutan <b>Organize favorites</b> (Organisera favoriter). Se 6.4 Start Favorite (Starta Favorit) för ytterligare information.
Kryssrutan Run this method now (Kör denna metod nu)	Markera den här kryssrutan om du vill starta metoden omedelbart efter att du har avslutat den här guiden.

Det ändrade dokumentet kan sparas under sitt nuvarande filnamn eller under ett annat namn.

Klicka på Save&Finish (Spara och slutför) för att spara metoden och stänga guiden Create/Edit a Method (Skapa/redigera en metod)



Klicka på den här knappen för att öppna dialogrutan Knappen Audit trail... Audit trail... där användaren, datumet, den utförda åtgärden och en kommentar visas. En förhandsgranskning av arbetsytan, metoden och prov-ID-listan för den aktuella och alla tidigare versioner kan visas genom att klicka på lämpliga knappar. Dessutom är det möjligt att spara en tidigare version av metoden under ett annat namn. Audit Trail User Date Action Comment 1 Administrator (Admin) 2 Administrator (Admin) 18.10.2005 15:57:39 Created (.wsp Modified) 18.10.2005 16:29:24 Modified (.wsp Modified) demo < Print preview Workspace... Method... Sample ID list OK Cancel Help Knappen Klicka på den här knappen för att öppna dialogrutan Signature... Signature... (Underskrift) där användare, datum, åtgärd (Underskrift) och kommentar visas. Den här knappen är endast tillgänglig när filen har signerats med hjälp av guiden Attach Signature (Bifoga underskrift). Se kapitel 8 Guiden Bifoga underskrift för ytterligare detaljer. Fältet När du anger ett lösenord för en metod kommer Method password metoden att vara lösenordsskyddad och kan inte ändras (Lösenord för metod) förrän rätt lösenord har angetts. Se detaljerad beskrivning nedan.

Ytterligare knappar i magellan Tracker:

# 4.4.1 Lösenordsskydd av metoder

#### Skydda metod med ett lösenord

Användare som har rättigheten att skapa och redigera metoder kan skydda metoder genom att tilldela dem ett lösenord.

Som standard har Administrators (Administratörer) och Application Specialists (Programspecialister) de nödvändiga rättigheterna.

När en ny metod skapas kan lösenordsskyddet ställas in på fliken **Save as** (Spara som) i guiden **Create/edit a method (Skapa en metod)**. Ange metodens namn, skriv lösenordet i redigeringsfältet **Method password (Lösenord för metod)** i (dialogrutans nedre högra hörn) och klicka på **Finish (Slutför)**.

Bekräfta lösenordet i dialogrutan **Method password (Lösenord för metod)** som visas och klicka på **OK**. Metoden sparas med lösenordsskydd.

Redan sparade, men inte lösenordsskyddade metoder kan skyddas genom att öppna metoden och ange lösenordet i fliken **Save as tab (Spara som)** enligt beskrivningen ovan.

#### Redigera lösenordsskyddade metoder

Om du vill öppna en lösenordsskyddad metod anger du det korrekta lösenordet för metoden. Metoden öppnas i **Edit mode (Redigeringsläge)**. Det är möjligt att redigera och spara metoden.

Om **Read only (Skrivskyddad)** väljs, öppnas metoden i skrivskyddat läge. Det är fortfarande möjligt att redigera metoden, men det är inte möjligt att spara ändringarna.

#### Ändra lösenordet på en lösenordsskyddad metod

För att ändra lösenordet för en lösenordsskyddad metod öppnar du metoden i Edit mode (Redigeringsläge) (lösenord krävs) och ange ett nytt lösenord i redigeringsfältet Method password (Lösenord för metod) i Fliken Spara som (Spara som). Klicka på Finish (Slutför).

Bekräfta lösenordet i dialogrutan **Method password ((Lösenord för metod)** och klicka på **OK**.

# 4.5 Metoder med flera plattor

Vid anslutning till ett INFINITE F50-instrument kan metoder med flera plattor definieras.

Antalet plattor som ska mätas kan ställas in i redigeraren för mätparametrar under metoddefinitionen.

deauwersert parameters - infante FSD		×
🔶 Lab Ware 🏦	🔷 🕈 Rate	Selecton
🔶 Flate	Pete definition: [GPE2004_Greener 56 Flat Transport.	Nothing selected
• Part of Plate	Une a start of the plane	
V Heasurements 2	* PatalPlan	
4 Absorbance		-
Actions E	A B B B B B B B B B B B B B B B B B B B	
N Station		
R, Nove Plate		
Kawtic 2	· 00000000000	
G Kinetic Cycle	" ••••••••••••••••••••••••••••••••••••	
A Kinetic Condition		8
Macelaeeus 2	↓ ▼ Absolutes	
Convert	Westmany Difference and Difference a	
<ul> <li>Visit (Timer)</li> </ul>	Viewerson Rom v	
incubation		
		~
Number of plates 3 2		
Help		
Carcel «««liack	CHOOSE MEASUREMENT	PARAMETERS
8 1990 Turns		

När två eller flera plattor ska mätas finns en räknare tillgänglig för att växla mellan plattorna i fönstret **Define Evaluation (Definiera utvärdering)** också.



#### 4. Create/Edit a Method Wizard (Guiden Skapa/redigera en metod)

No.         No. <th>941,81 17 941,82 17 941,82 941,83</th>	941,81 17 941,82 17 941,82 941,83
ST(1)         SM(2)         SM(2) <th< td=""><td>SW1,02 V1 SW1,02</td></th<>	SW1,02 V1 SW1,02
STE2         SME, 20         S	501,03
*** 515 51 51 51 51 51 51 51 51 51 51 51 51	In part 1
515.4 SM1_5 SM1_12 SM1_21 SM1_21 SM1_21 SM1_21 SM1_22 SM1_24 SM1_20 SM1_24 SM1_20 SM1_24 SM1_20 SM1_24 SM1_	50m
	Raphann 1111 B. Tanacter
ST1_3         SM(-6)         SM(-3)         SM(-2)         SM(-3)         SM(-3)         SM(-4)         SM(-3)         SM(-3)         SM(-4)         SM(-4)         SM(-3)         SM(-3)         SM(-4)         SM(-4) </td <td>CM Internet</td>	CM Internet
PCT         SML_2         SML_35         SML_25         SML_35	50/1_87 1/1
HOT 397, B 397,	5841_68 1/1

Alla andra parametrar kan definieras på samma sätt som för metoder med en enstaka platta. När du startar en metod med flera plattor visas följande dialogruta:

Run measurement		
Measurement of	plate 🚺 of 3	
Sample IDs list:		Insert
	STOP PAUSE RUN	
		Plate in

Sample IDs List (Prov-ID-lista)	För varje enstaka platta kan en prov-ID-lista väljas.
Knappen STOP (Stopp)	Mätningen stoppas. Data för de redan uppmätta plattorna visas i det skapade arbetsytan och det är inte möjligt att lägga till mätningar av ytterligare plattor.
Knappen PAUSE (Paus)	Mätningen pausas. Data för de redan uppmätta plattorna visas i det skapade arbetsytan, men det är möjligt att lägga till mätningar av ytterligare plattor.
Knappen RUN (Kör)	Mätning av nästa platta påbörjas.
Plate in/ Plate out (Platta in/platta ut)	Plattan flyttas in i eller ut ur instrumentet.

När du kör en metod med flera plattor kan du pausa mätningen mellan två plattor, spara de uppnådda resultaten och fortsätta mätningen för de återstående plattorna vid ett senare tillfälle. Om **Automatic save workspace (Spara arbetsyta automatiskt)** väljs, sparas arbetsytan efter varje enstaka mätning.

# 5. Create/Edit a Sample ID List Wizard (Guiden Skapa/redigera en prov-ID-lista)

# 5.1 Inledning

Create/edit a sample ID list wizard (Guiden Skapa/redigera en prov-ID-lista) används för att skapa nya provlistor eller redigera befintliga provlistor.

# Sammanfattning av arbetsflöde

Starta guiden **Create/edit a sample ID list (Skapa/redigera en prov-ID-lista)**. Efter en välkomstsida visas sidan **File selection (Filval)**. Välj ett av alternativen: **Create new (Skapa ny)** prov-ID-lista eller **Edit (Redigera)** en befintlig prov-IDlista och spara ID-listan med hjälp av sidan **Save as (Spara som)**.

# 5.2 Create/Edit a Sample ID List (Skapa/redigera en prov-ID-lista)

Alternativknappen Create new (Skapa ny)	En ny prov-ID-lista kan skapas.
Alternativknappen <b>Edit</b> (Redigera)	En befintlig prov-ID-lista kan ändras.
Kombinationsrutan <b>Show</b> (Visa)	<ul> <li>Filerna visas i enlighet med det valda urvalet:</li> <li>All files (Alla filer)</li> <li>My files (Mina filer): Det här alternativet är tillgängligt om användaradministrationen är aktiverad (alltid aktiverad i magellan Tracker)</li> </ul>
Listan <b>Filename</b> (Filnamn)	Den prov-ID-lista som ska redigeras måste väljas. Fältet <b>Remarks</b> (Anmärkningar) bredvid varje filnamn innehåller - om det anges - en kort beskrivning av prov- ID-listan. Alla prov-ID-listor som finns tillgängliga i standardkatalogen för prov-ID-listor och som uppfyller de filterkriterier som valts i kombinationsrutan show visas.
Knappen <b>Print Preview</b> (Förhandsgranskning av utskrift)	En förhandsgranskning av den valda prov-ID-listan visas och en utskrift kan påbörjas.

File Selection (Filval) innehåller följande element:

# 5.2.1 Skapa ny prov-ID-lista)

I fönstret **Select Plate Type (Välj plattyp)** kan antalet prov-ID:n per brunn och den plattypen som ska användas väljas. Alternativet **Import from external sample ID list (Importera från extern prov-ID-lista)** möjliggör sökning efter olika typer av prov-ID-filer och importerar dem till **magellan**.



Anmärkning Vissa prov-ID-listor kan innehålla ytterligare information, t.ex. koncentrationer och spädningar, som kan påverka utvärderingen av mätdata.



Select plate type Plate type Plate type Plate type from method Method1.mth Load Standard plate types S6 well plate Plate type from pdf-file CL196fb.pdf Import Import	Settings Sample ID's per well ② @ne ③ Iwo ③ Ihree Printout ④ horiz. Table ④ vert. Table ④ vert. Table ④ Matrix
Help Cancel <<< Back	ENTER SAMPLE IDs

Dialogrutan Select plate type (Välj plattyp) innehåller följande element:

Settings (Inställningar)	
Alternativknapparna <b>Sample IDs per well</b> (Prov-ID:n per brunn)	Ange det antal prov-ID:n som krävs per brunn (maximalt tre kan anges).
Alternativknapparna Printout (Utskrift)	Definiera layouten för utskrift av prov-ID-listan.
Alternativknappen <b>Plate type</b> (Plattyp)	Fyra alternativknappar definierar vilken plattyp som ska väljas:
	<ul> <li>Plate type from method (Plattyp från metod) Välj en metod för att läsa in plattlayouten. Klicka på Load (Ladda) för att söka efter befintliga metodfiler. Dialogrutan Windows<sup>TM</sup> Open (Öppna) visas. Välj önskad fil och klicka på Open (Öppna).</li> <li>Standard plate type (Standard plattyp) Plattstyper med standardgeometri kan väljas för att definiera plattformatet. När den är aktiverad kan en standardgeometrisk plattyp väljas från en tillhörande rullgardinsmeny.</li> </ul>
	<ul> <li>Plate type from pdf-file (Plattyp från PDF-fil) En fördefinierad pdf-fil (SUNRISE-instrument) eller en pdfx-fil (INFINITE F50-instrument) kan användas för att definiera plattformatet.</li> </ul>
	<ul> <li>Import from external Sample ID list (Importera från extern prov-ID-lista)</li> <li>En extern filtyp kan väljas för import som prov-ID-lista.</li> <li>Se kapitel 5.2.3 Importera en prov-ID-lista.</li> </ul>

Knappen **Enter sample IDs (Ange prov-ID:n)** är endast aktiverad om de nödvändiga inställningarna har definierats (metod har laddats, standardplattyp eller PDF-fil har valts).

# 5.2.2 Import/Edit a Sample ID List (Importera/redigera en prov-ID-lista)

Efter att ha valt plattyp, visas fönstret **Import/edit a sample ID list** (Importera/redigera en prov-ID-lista):

- Verktygsfält med knapparna Settings (Inställningar), Edit (Redigera), Autofill (Autofyll), Delete (Radera), Undo/Redo (Ångra/Gör om), Print format (Skriv ut format) Print preview (Förhandsgranskning av utskrift), och Zooming (Zooma).
- Plattvy med en schematisk layout av en mikroplatta.
- Prov-ID-lista (när Plate type from method (Plattyp från metod) har valts)
- Dialogruta för att infoga prov-ID:n (när Standard plate type (Standard plattyp) har valts)

Om en ny prov-ID-lista ska skapas tilldelas inga prov-ID:n till brunnarna. En brunn kan väljas genom att klicka på brunnen.

Alla nödvändiga prov-ID:n måste anges antingen manuellt eller med hjälp av knapparna i verktygsfältet. De kan även läsas av den handhållna streckkodsläsaren för prov-ID:n. Prov-ID:n kan definieras och redigeras i den definierade mikroplattstypen. Upp till tre prov-ID:n kan anges i varje brunn. Prov-ID:n kan flyttas eller kopieras.

När en metod har valts visas listan Lista över alla prover från plattans layout (Lista över alla prover från plattlayouten) på vänster sida. Endast prover utan replikat visas i listan. Prov-ID:n kan anges antingen i listan eller på plattan. Alla ändringar visas omedelbart.

Knapr	bar i	verktvo	asfältet

Knappen Settings (Inställningar)	Knappen öppnar <b>Sample ID List Settings (Inställningar för prov-ID-listan)</b> där antalet prov-ID:n per brunn kan väljas.
Knappen Edit (Redigera)	Knappen <b>Edit (Redigera)</b> öppnar dialogrutan <b>Edit Well</b> (Redigera brunn). Prov-ID:n för den markerade brunnen kan redigeras. Se nedan för mer information.
Knappen Autofill (Autofyll)	Hela plattan eller en del av plattan kan väljas för att automatiskt fyllas med prov-ID:n. När du har valt motsvarande brunnar klickar du på <b>Autofill (Autofyll)</b> för att öppna dialogrutan <b>Autofill Selection (Autofyll urval)</b> . Se nedan för mer information.
Knappen Delete (Radera)	Hela plattan eller en del av plattan kan väljas. Den här funktionen raderar prov-ID:n för de valda brunnarna omedelbart. Ingen dialogruta begär en bekräftelse. Klicka på <b>Delete (Radera)</b> eller tryck på <b>DEL</b> för att radera prov-ID:n i de valda brunnarna.
Knappen Undo (Ångra)	Ångrar den senaste åtgärden. Klicka på <b>Undo (Ångra)</b> eller tryck på <b>CTRL-Z</b> för att ångra en tidigare åtgärd.
Knappen Redo (Gör om)	Upprepar den åtgärd som utfördes före ångra-steget. Klicka på <b>Redo (Gör om)</b> eller tryck på <b>CTRL+Y</b> för att göra om en åtgärd som ångrades.

Print format (Skriv ut format)	<ul> <li>Välj layout för utskrift av prov-ID-listan:</li> <li>horiz. Table (Horisontell tabell)</li> <li>vert. Table (Vertikal tabell)</li> <li>Matrix (Matris)</li> </ul>
Print Preview… (Förhandsgransknin g av utskrift)	Öppnar förhandsgranskningen av den utskrivna prov-ID- listan.
Knappen 10 %	Detta ställer in plattlayoutvyn på 10 %.
Knappen 100 %	Detta ställer in plattlayoutvyn på 100 %.
Kryssrutan Zoom	Om brunnsplattan består av ett stort antal brunnar kan de individuella brunnstilldelningarna inte visas i 100 %- zoomläget. I ett sådant läge tillåter knappen <b>Zoom mode</b> ( <b>Zoom-läge</b> ) att det markerade området zoomas in. Om den är aktiverad kan användaren välja ett område att zooma in i genom att markera det önskade layoutområdet. Klicka på höger musknapp för att zooma ut till 100 %.



# Sammanhangsberoende meny för plattlayouten

Genom att högerklicka på en brunn i plattlayouten visas en sammanhangsberoende meny för de markerade brunnarna. Följande kommandon är tillgängliga:

Meny	Beskrivning
Summary (Sammanfattning)	Dialogrutan <b>Summary (Sammanfattning)</b> visas. Se nedan för mer information.
Cut (Klipp ut)	Prov-ID:n för de markerade brunnarna tas bort från brunnarna och kopieras till urklippet.
Copy (Kopiera)	Prov-ID:n för de markerade brunnarna kopieras till urklippet.
Paste (Klistra in)	Prov-ID:n som kopierats till urklippet klistras in i de markerade brunnarna.
Delete (Radera)	Klicka på <b>Delete (Radera)</b> eller tryck på <b>DEL</b> för att radera prov-ID:n för de markerade brunnarna.
Copy in ASCII- Format (Kopiera i ASCII-format)	Prov-ID:n för de markerade brunnarna kopieras till urklippet i ASCII-filformat. På så sätt kan de enkelt klistras in i andra program (t.ex. Excel, Notepad).
Paste in ASCII- Format (Klistra in i ASCII-format)	Prov-ID:n kan klistras in från urklippet i ASCII-filformat från andra program. Se nedan för mer information.
Redigera	Alternativet <b>Edit (Redigera)</b> öppnar dialogrutan <b>Edit Well</b> (Redigera brunn). Prov-ID för den markerade brunnen kan redigeras. Se nedan för mer information.
Autofill Selection (Autofyll urval)	Hela plattan eller en del av plattan kan väljas för att automatiskt fyllas med prov-ID:n. Efter att ha valt motsvarande brunnar väljer du alternativet Autofill (Autofyll) och dialogrutan Autofill Selection (Autofyll urval) visas. Se nedan för mer information.



Anmärkning När du redigerar en prov-ID-lista kan du använda funktionen dra-ochsläpp. Använd vänster musknapp för att välja och dra elementet.

#### Edit well (Redigera brunn)...

Dialogrutan **Edit Well (Redigera brunn)** kan nås via den sammanhangberoende menyn eller knappen i verktygsfältet.

Prov-ID:n för alla brunnar på en mikroplatta kan definieras. Maximalt tre prov-ID:n per brunn är möjliga.

De önskade prov-ID:n för den markerade brunnen måste anges i dialogrutan **Edit Well (Redigera brunn)**. Detta måste upprepas för varje brunn som ska fyllas med prov-ID:n. Navigationsfunktionerna i dialogrutan **Edit Well (Redigera brunn)** kan användas för att underlätta arbetet. Klicka på **OK** för att stänga dialogrutan, och prov-ID:n för den senast markerade brunnen sparas.

Dialogrutan Edit Well (Redigera brunn) innehåller följande element:

Knapparna Left (Vänster), Right (Höger), Up (Upp) och Down (Ner)	Riktningsknapparna kan användas för att välja en annan brunn på plattan samtidigt som dialogrutan förblir öppen och ändringarna i den aktuella brunnen sparas. <b>Piltangenterna</b> kan också användas för denna typ av navigering.
Kryssrutan On Enter move in vertical direction (Vid Enter förflytta i vertikal riktning)	När du klickar på <b>Enter</b> eller <b>Tab</b> flyttas inmatningsfältet till nästa brunn. Använd kryssrutan för att ange rörelseriktningen, vertikalt eller horisontellt över plattan.
Textfältet Edit sample IDs (Redigera prov- ID:n)	I varje brunn kan upp till tre prov-ID:n anges. Listan innehåller därför tre <b>Sample ID (Prov-ID)</b> -textfält och kryssrutan <b>Autoincrement (Automatisk ökning)</b> som är associerad med vart och ett av dem. Ett prov-ID för den markerade brunnen kan anges i textfälten. Prov-ID kan innehålla bokstäver, siffror eller en kombination av båda. Till exempel: Smp 1. Upp till 100 alfanumeriska tecken kan anges. Hela plattan kan fyllas med prov-ID.
Kryssrutan Autoincrement (Automatisk ökning)	Kryssrutan Autoincrement (Automatisk ökning) kan markeras om brunnarna för prov-ID:n ska numreras automatiskt. Till exempel: Om startvärdet är Smp1 kallas nästa brunn automatiskt för Smp2, med hjälp av Enter/Tab eller pilknapparna. Med hjälp av Autoincrement (Automatisk ökning) kan prov-ID:n tilldelas till brunnarna på ett numeriskt sammanhängande sätt. Ange bara det ursprungliga värdet i den första brunnen. Det angivna ID:t kan också vara alfanumeriskt, även om det måste sluta med ett nummer för att möjliggöra t.ex. sammanhängande numrering av ytterligare brunnar: Om John1 anges i den första brunnen kommer John2 automatiskt att visas när du flyttar till den andra brunnen.

### Autofill Selection (Autofyll urval)

Dialogrutan Autofill Selection (Autofyll urval) kan nås via den sammanhangberoende menyn eller knappen i verktygsfältet.

Denna dialogruta gör det möjligt för användaren att automatiskt fylla de valda brunnarna med prov-ID:n.

Välj hela plattan eller delar av plattan och ange önskade prov-ID:n och ökningsinställningar för de markerade brunnarna i dialogrutan Autofill Selection (Autofyll urval). Klicka på OK för att stänga dialogrutan och för att fylla de valda brunnarna med de definierade prov-ID:n. I varje brunn kan upp till tre prov-ID:n anges. Listan innehåller därför tre textfält för **Sample ID** (prov-ID) och en kryssruta för **Autoincrement** (Automatisk ökning), ett textfält för **Startindex** och en kryssruta för **Vertical** (Vertikal) som är associerad med vart och ett av dem.

Dialogrutan Autofill Selection (Autofyllt urval) innehåller följande element:

Kryssrutan <b>Use ID</b> (Använd ID)	Kryssrutan <b>Use ID</b> (Använd ID kan endast användas om ett identiskt antal prov-ID:n har valts. Krysset avgör om funktionen för automatisk ökning tillämpas på det valda prov-ID:t.
Textfälten <b>Sample ID</b> (Prov-ID)	Prov-ID:n för den markerade brunnen kan anges i textfältet <b>Sample ID</b> (Prov-ID). Prov-ID:n kan innehålla bokstäver, siffror eller en kombination av bokstäver och siffror.
Kryssrutan <b>Autoincrement</b> (Automatisk ökning)	Kryssrutan <b>Autoincrement</b> (Automatisk ökning) måste vara markerad om prov-ID:n för brunnarna ska numreras automatiskt (ökas med 1).
Textfältet Startindex	Textfältet <b>Startindex</b> är endast redigerbart om <b>Autoincrement</b> (Automatisk ökning) är vald. Startvärdet för den automatiska ökningen av prov-ID måste anges.
Kryssrutan <b>Vertical</b> (Vertikal)	Kryssrutan <b>Vertical</b> (Vertikal) kan endast användas om kryssrutan <b>Autoincrement</b> (Automatisk ökning) är markerad. Om den här kryssrutan är markerad kommer prov-ID:n att anges och numreras i vertikal riktning. Om den är avmarkerad kommer prov-ID:n att anges och numreras i horisontell riktning.

#### Well Summary (Sammanfattning av brunn)

Dialogrutan **Well Summary (Sammanfattning av brunn)** kan nås via den sammanhangsberoende menyn.

Den här dialogrutan ger en översikt över alla definierade prov-ID:n och de inställningar som är relaterade till prov-ID-listan för en vald brunn.

Dialogrutan Summary... (Sammanfattning) innehåller följande element:

Information tree (Informationsträd)	Informationsträdet ger en översikt över alla definierade brunnsbaserade parametrar för den valda brunnen. Trädet fylls med den information som finns tillgänglig. För prov-ID-listor visas prov-ID:n, pipetteringsstatus osv.
Knapparna Left (Vänster), Right (Höger), Up (Upp) och Down (Ner)	Riktningsknapparna kan användas för att välja en annan brunn på plattan samtidigt som dialogrutan förblir öppen.
Expand All/ Shrink All (Expandera allt/förminska allt)	Den här knappen visar informationsträdet expanderat till den högsta nivån eller förminskat till den första nivån.

#### Paste in ASCII-Format (Klistra in i ASCII-format)

Funktionen **Paste in ASCII-Format (Klistra in i ASCII-format)** finns tillgänglig via den sammanhangsberoende menyn.

Den här funktionen klistrar in innehållet i utvalda data i ASCII-format från urklippet till **magellan**.

Funktionen gör det möjligt att överföra data från andra programvaror (t.ex: Excel, Notepad) till **magellans** plattlayout.

Först måste de relevanta data väljas i det andra programmet och kopieras till urklippet. Raderna måste skiljas åt med radmatning och kolumnerna med tabbstoppar. Detta görs automatiskt i Excel när du markerar flera celler.

Funktionen **Paste in ASCII-Format (Klistra in i ASCII-format)** klistrar in de kopierade data ASCII-format från urklippet till brunnarna. Data infogas alltid från position A1.

# 5.2.3 Importera en prov-ID-lista

Prov-ID-listor som har skapats med andra program eller prov-ID-listor som har definierats med andra former av analysutrustning måste importeras till **magellan** innan de används.



#### FÖRSIKTIGHET VISSA PROV-ID-LISTOR KAN INNEHÅLLA YTTERLIGARE INFORMATION, T.EX. KONCENTRATIONER OCH SPÄDNINGAR, SOM KAN PÅVERKA UTVÄRDERINGEN AV MÄTDATA.

Klicka på knappen **Import (Importera)** i dialogrutan **Sample ID list settings (Inställningar för prov-ID-lista)** för att söka efter befintliga externa prov-ID-filer. Dialogrutan Windows<sup>TM</sup> Open (Öppna) visas.

I fönstret visas en lista över File names (Filnamn) som matchar det valda File type (Filtypen). File type (Filtypen) kan väljas från en tillhörande rullgardinsmeny.

Följande externa filtyper stöds:

- Easy-Files .esy
- Tecan-files .tpl
- DD1-filer .dd1
- · Hamilton-Files .pro
- APL-filer .apl
- Gemini-filer .csv
- Anpassade formatfiler .txt

Välj den begärda filen som ska importeras och klicka på knappen **Open (Öppna)**. Dialogrutan Windows<sup>™</sup> **Open (Öppna)** stängs och filen laddas.

Bekräftelse av dialogrutan *Sample ID List Setting* (Inställning av prov-ID-lista) med Next (Nästa) kommer att öppna fönstret *Import/Edit a Sample ID List* (Importera/redigera en prov-ID-lista) med prov-ID:n som visas.

Knappen Import (Importera) kan också nås via dialogrutan Insert Sample ID List (Infoga prov-ID-lista) som finns i fönstret Start Measurement (Starta mätning) (Knappen Insert (Infoga)) och plattvyn Utvärdera resultat (Edit (Redigera) i verktygsfältet).



# FÖRSIKTIGHET

KONTROLLERA ATT RÄTT ID-LISTA HAR VALTS, ANNARS KOMMER RESULTATEN ATT TILLDELAS EN ANNAN LISTA.





Anmärkning När du importerar prov-ID-listor i .tpl, .dd1 eller anpassade format kan ytterligare information som ändrar utvärderingsdefinitionerna

inkluderas.

#### Anmärkning

När du importerar prov-ID-listor med spädningar får spädningen inte vara större än 1:1e+12 (magellan Standard och magellan Tracker).



#### Easy-Files .esy

Dessa filer består av rader som innehåller prov-ID och brunnsnamn.

Well name (Brunnens namn):	alfanumerisk, max. 3 tecken
Sample-ID (Prov-ID):	alfanumerisk, max. 20 tecken

Raderna skiljs åt med CR + LF.

#### Exempel

C01

B01

A01 0001

H02 0002

G02 0003

F02 0004

E02 0005

#### Tolkning

När magellan behandlar en .esy-fil

- Antalet prov-ID:n är satt till 1
- Prov-ID anges som prov-ID i brunnen

#### **Tecan-files** .tpl

Dessa filer (<PLATEID>.TPL) består av rader som innehåller parametrar separerade med semikolon. Formatet är:

H;date;time (H;datum;tid)

D;testno;sample-ID + control-name;position;pre-dilution;errors (D;testnr;prov-ID + kontrollnamn;position;före spädning;fel)

L;

Filen börjar med ett H och slutar med ett L. Fälten definieras på följande sätt:

date;time;testno;sample-ID (datum, tid, testnr, prov-ID)	samma som i jobblistan (BES3)
control name (kontrollnamn)	alfanumerisk, max 8 tecken Anmärkning! Denna inmatning innehåller prov-ID eller namnet på en kontroll
position	alfanumerisk, max 3 tecken
pre-dilution (för-spädning)	numerisk, max. 8 tecken plus 1 decimal (separerade med ett ".")

errors (fel)	alfanumerisk, max 3 tecken Om ett fel inträffade under pipetteringen visas ett av följande tecken här:		
	Första teckn	et	
	<mellanslag:< td=""><td>&gt;</td><td>ingen vätska</td></mellanslag:<>	>	ingen vätska
	L:	ingen	eller för lite vätska
	Andra teckne	ət	
	<mellanslag:< td=""><td>&gt;</td><td>ingen streckkod</td></mellanslag:<>	>	ingen streckkod
	M:	manu	ellt angiven streckkod
	Tredje teckn	et	
	<mellanslag< td=""><td>&gt;</td><td>ingen överskridande av timer</td></mellanslag<>	>	ingen överskridande av timer
	O:	timer	överskridande

Raderna skiljs åt med CR + LF.

#### Exempel

H;29-02-94;12:47:11

D;1;NEG;A1;1.0;

D;1;POS;B1;1.0;

D;1;0001;C1;1.0

L;

#### Tolkning

När magellan behandlar en .tpl-fil

- Antalet prov-ID:n är satt till 1
- Prov-ID eller kontrollnamn är inställt som prov-ID i brunnspositionen
- Felmarkeringar lagras för varje brunn
- För-spädning används som spädning av brunnspositionen

#### DD1-filer .dd1

Dessa filer består av rader som innehåller parametrar med en fast längd. Formatet är:

Första raden: Datum (DD-MM-YY)

Andra raden: Tid (HH:MM:SS)

Fält	Pos	Längd	Format/Beskrivning
TestNr	1–3	3	nummer (001–049)
TestName	4–23	20	Alfanumerisk
Prov-ID	24–35	12	Alfanumerisk
ControlName	36–43	8	Alfanumerisk
PlateID	44–51	8	Alfanumerisk
PlatePos	52–55	4	Alfanumerisk
Fel	56–61	6	Alfanumerisk

Raderna skiljs åt med CR + LF.

# Exempel

13-04-1999 11:01:32			
953test1 KBR		KBR3	H1
953test1 KBR		KBR3	G1
953test1 KBR		KBR3	C1
953test1 KBR		KBR3	B1
953test1 KBR		KBR3	A1
953test1 KBR	0001	KBR3	H2
953test1 KBR	0001	KBR3	G2
953test1 KBR	0001	KBR3	F2
953test1 KBR	0001	KBR3	E2
953test1 KBR	0001	KBR3	D2

#### Tolkning

När magellan behandlar en .dd1-fil

- Antalet prov-ID:n är satt till 1
- <Testname> anger magellan metodnamnet i brunnen
- <PlatePos> anger brunnen
- <SampleID> eller <ControlName> anger magellan prov-ID.
- <Errors> anger pipetteringsflaggan: L Vätskefel, C: Koagelfel, E: manuellt inmatat

När en .dd1-fil läses kommer den aktuella layouten att ändras med hjälp av den layout som definieras i de olika metoderna där namnen anges i .dd1-filen.

#### Hamilton-Files .pro

Dessa filer består av rader som innehåller prov-ID och brunnens namn.

Det finns också information om sidhuvudet.

Well name (Brunnens namn): alfanumerisk, max. 3 tecken

Sample-ID (Prov-ID): alfanumerisk, max. 20 tecken

Raderna skiljs åt med CR + LF.

### Exempel

14ESP-AK-Bommeli	-> Sidhuvud
08403009070	-> Sidhuvud
SendInfoHost=00	-> Sidhuvud
SendInfoPrnt=00	-> Sidhuvud
SendInfoDisk=00	-> Sidhuvud
A0109111111110	

A02090325228840

A03090325228930

A04090325235730

A05090336975200



#### Tolkning

När magellan behandlar en .pro-fil

- Antalet prov-ID:n är satt till 1
- Prov-ID är inställt som prov-ID i brunnen
- Information om sidhuvud ignoreras

#### **APL-filer** .apl

Beskrivning av .apl-filer skapade av MiniSwift 1.3 eller senare, eller Minilyser 5.30 eller senare. Filformatet definieras på följande sätt:

Fält	Pos	Längd	Format/Beskrivning
PlateID	1–8	8	Alfanumerisk
PlatePos	9–12	4	Alfanumerisk
TestNumber	13–15	3	Antal
TestName	16–27	12	Alfanumerisk
SampleID	28–39	12	Alfanumerisk
ControlName	40–47	8	Alfanumerisk
Result_1	48–55	8	Alfanumerisk
Result_2	56–63	8	Alfanumerisk
Errors	64–69	6	Alfanumerisk
Date	70–79	10	Alfanumerisk
Time	80–87	8	Alfanumerisk
CutOff	88–95	8	Alfanumerisk

Raderna skiljs åt med CR + LF.

#### Exempel

990520S1A1	2	test1_2	BLANK		05-20-199919:12:53
990520S1B1	2	test1_2	BLANK		05-20-199919:12:53
990520S1C1	2	test1_2	NC1	LR	05-20-199919:12:53
990520S1D1	2	test1_2	PC1		05-20-199919:12:53
990520S1E1	2	test1_2	0000000010		05-20-199919:12:53
990520S1F1	2	test1_2	0000000010	V	05-20-199919:12:53

#### Tolkning

När magellan behandlar en .apl-fil

- Antalet prov-ID:n är satt till 3
- <Testname> anger magellan metodnamnet i brunnen
- <PlatePos> anger brunnen
- <SampleID> anger den magellan Prov-ID 1
- <ControlName> anger den magellan Prov-ID 2
- <Errors> anger den magellan Prov-ID 3
- <Errors> anger pipetteringsflaggan: pipetteringsfel L, B, N i position 1 är okritiska för prover men kritiska för kontroller, fel i verifieringsläsning och streckkod är också okritiska.



#### Gemini-filer .csv

Beskrivning av .csv-filer som skapats av Gemini 3.40 och 3.60. Följande är ett exempel på en utdatafil i Excel-format:

Position	Tube-ID	ScanError	SRCRack	SRCPos	SRCTubeID	Volym
Error	SRCRackID	GridPos	SiteOnGrid	TipNumber	DetectVol	Time
	Version 2.00	3091	Microplate	12	8	
0	9900001	18	1			3.8.1998
1		0	Tube 13*85mm	1	0001	10
0	013_000064	1	1	1		11:55:42
3		0	Tube 13*85mm	2	0002	20
0	013_000064	1	1	2		11:55:42
5		0	Tube 13*85mm	3	0003	30
0	013_000064	1	1	3		11:55:42
7		0	Tube 13*85mm	4	0004	40
0	013_000064	1	1	4		11:55:42

Filformatet definieras på följande sätt:

Den första raden innehåller kolumnrubrikerna, som är samma för alla format, utom för vissa format som endast stöder versaler. Den andra raden visar information om det stället/stripsstället som filen skapades för. Endast kolumnerna Tube ID, SRCRack, SRCPos, SRCTubeID, Error, SRCRackID, GridPos, SiteOnGrid och Time är giltiga.

Med Gemini 3.60 innehåller SRCPos antalet x brunnar, SRCTubeID antalet y brunnar och Volume innehåller den rad som ska läsas.

Den tredje och följande raderna är poster för de använda brunnarna i ett ställ/de använda rören i ett stripsställ.

Kolumnerna innehåller följande information (Column\Range\Explanation):

- Position\1 .. antal brunnar\Position för brunnen i stället. Numreringssekvensen är bak till fram och sedan från vänster till höger. (1 = A1, 2 = B1, ...) Om mikroplattan har pipetterats i stående läge (t.ex. dimensioner 8x12, iso. 12x8) är numreringssekvensen vänster till höger och sedan fram till bak. (1 = H1, 2 = H2, ...)
- TubeID\Max. ra32 tecken\Denna kolumn fylls endast i för stripsställ och när rörens streckkoder skannades med ett POSID-kommando. I rad 2 visar denna kolumn versionen av Gemini.
- ScanError\0.. 4294967295\Denna kolumn innehåller möjliga skanningsfel. Observera tabellen med felkoderna nedan. I rad 2 visar denna kolumn serienummer av Genesis.
- SRCRack\Max. 32 tecken\Här rapporterar Gemini konfigurationsnamnet för källstället för dispenseringskommandon. I rad 2 visar denna kolumn konfigurationsnamnet för det ställ för vilket utdatafilen skrevs.
- SRCPos\1 .. antal brunnar\Denna kolumn anger positionen i källstället, varifrån vätskan kom, för dispenseringskommandon.
- SRCTubeID\Max. 32 tecken\Om källan för ett dispenseringskommando var ett stripsställ rapporterar Gemini här streckkoden för källröret.



- Volym\-7158278 .. +7158278\Denna kolumn innehåller volymen i µl, som dispenserades (positivt värde) eller aspirerades (negativt värde).
- Error\0 .. 4294967295\Denna kolumn innehåller möjliga pipetteringsfel. Observera tabellen med felkoderna nedan. I rad 2 visar denna kolumn skanningsfelen.
- SRCRackID\max. 32 tecken\För dispenseringskommandon, är detta källställets streckkod. I rad 2 visar denna kolumn streckkoden för det ställ för vilket utdatafilen skrevs.
- GridPos\1..67 Detta är rutnätspositionen för källstället för dispenseringskommandon, .
- SiteOnGrid\1 ... 128\Här visar Gemini y-positionen (platsen på hållaren) på källstället för dispenseringskommandon.
- TipNumber\1 ... 8\Här rapporterar Gemini spetsnumret som användes för pipettering till/från denna brunn.
- DetectVol\0 .. +7158278\Om vätskedetektering användes, anges här den upptäckta volymen i brunnen före aspirering-/dispenseringskommandot.
- Tid\...\Detta är tiden (format hh:mm:ss) när brunnen användes.

Raderna skiljs åt med CR + LF. Kolumnerna skiljs åt med antingen kommatecken eller semikolon. Kolumnavskiljare identifieras från den första rubrikraden.

#### Tolkning

Antalet rader separerade med en ny rad räknas. Detta antal reducerat med 2 används för att definiera plattformatet, som kan vara 8x12, 16x24, 24x36, 32x48 (rader x kolumner).

Med Gemini 3.60 extraheras formatinformationen från den andra raden, SRCPos och SRCTubeID. Flera rader per brunn är möjliga. Siffran på den andra raden, Volume (Volym), anger vilken rad som ska användas. Om den angivna raden inte finns, kan prov-ID-listan inte läsas. Om någon rad innehåller ett felnummer som är större än noll, anges antingen ett kritiskt eller ett okritiskt fel, beroende på felnumret.

- Position: Well position (t.ex. 1 = A1, 2 = B1, ...)
- Tube-ID: ignorerad
- Skanningsfel: Om felnumret är större än noll anges ett kritiskt fel
- SRCRack: prov-ID 2
- SRCPos: om SRCTubeID är tomt, används som prov-ID 1 (SRCRackID + "\_" + SRCPos)
- SRCTubeID: om det inte är tomt används det som prov-ID 1
- Volume (Volym): ignorerad
- Error (Fel): prov-ID 3, om felnumret är större än noll anges ett kritiskt fel
- SRCRackID: om SRCTubeID är tomt, används som prov-ID 1 (SRCRackID + "\_" + SRCPos)
- GridPos: ignoreras
- TipNumber: ignoreras
- DetectVol: ignoreras
- Time: ignoreras



#### Anpassade formatfiler \*.\*

När du importerar filer i anpassade format visas dialogrutan **Custom Format** (Anpassat format). Dialogrutan **Custom Format (Anpassat format)** innehåller följande element:

Grupprutan <b>Plate type (Plattyp)</b>	I grupprutan Plate Type (Plattyp) kan plattformatet anges. Om formatet redan har angetts är redigeringsfälten inaktiverade.
Rader	Ange antalet rader.
Kolumner	Ange antalet kolumner.
Grupprutan File format (Filformat)	l grupprutan File format (Filformat) kan formatet för den angivna filen definieras.
File extension (Filtillägg)	I fältet <b>File extension (Filtillägg)</b> visas filtillägget och fältet är inaktiverat.



#### Anmärkning Om du använder den här funktionen i guiden Create/Edit a method (Skapa/redigera en metod) aktiveras fältet för filtillägg och filtillägget måste anges.

Kombinationsrutan	Välj den kolumnavskiljare som används i filen.
<b>Kolumnavskiljare</b>	Radavgränsare måste vara CR+LF.
Börja med raden	Ange ett tal större än 1 om rubrikrader ska ignoreras.

#### Exempel

Anta följande inställningar:

Plate Type (Plattyp):

- 2 rader
- 3 kolumner

#### Format

- kolumnavskiljare = Tabulator
- börja med raden = 2
- kolumn 1 = Plattposition (A1, A2, ...)
- kolumn 2 = prov-ID 1
- kolumn 3 = pipetteringsstatus

# Exempel:

Exempel:		
**Tecan anpassad fil**		
A1	BLANK	0
A2	NC	0
A3	SM001	0
B1	SM002	3
B2	SM003	0
B3	SM004	0



# 5.2.4 Spara prov-ID-listan

Klicka på Next (Nästa) i fönstret Import/Edit a Sample ID List (Importera/redigera en prov-ID-lista) och fönstret Save as (Spara som) öppnas, där den skapade eller ändrade prov-ID-listan kan sparas:

Fältet <b>Filename</b> (Filnamn)	Det här fältet är avsett för att ange eller ändra filnamnet för prov-ID-listan (.smp).
Fältet File remarks (Filkommentarer)	Fältet <b>File remarks</b> (Filkommentarer) gör det möjligt att lägga till några kommentarer till filen, som kommer att visas i till exempel dialogrutorna <b>Open file</b> (Öppna fil).
Fältet <b>Audit trail</b> comment (Audit trail kommentar)	Fältet <b>Audit trail comment</b> (Audit trail kommentar) gör det möjligt att lägga till några kommentarer, som lagras i Audit trail (endast tillgängligt för <b>magellan</b> Tracker).
Knappen Audit trail	Visar audit trail för prov-ID-listan (endast tillgänglig i <b>magellan</b> Tracker). Se 4.4 Spara metoden - Audit trail för mer information.

Knappen Save&Finish (Spara och slutför) är endast aktiverad när ett dokument har ändrats. Det ändrade dokumentet kan sparas under sitt nuvarande filnamn eller under ett annat namn.

Klicka på **Save&Finish** (Spara och slutför) för att spara prov-ID-listan och stänga guiden **Create/Edit a Sample ID List** (Skapa/redigera en prov-ID-lista).
# 6. Start Measurement Wizard (Guiden starta mätning)

# 6.1 Inledning

#### I fönstret Wizard List (Guide-lista) klickar du på Start Measurement (Starta

mätning) för att öppna guiden. Klicka på Continue (Fortsätt) bå välkomstsidan och dialogrutan Select a file (Välj en fil) visas:

Select a File		×	
Obtain Raw Data			
Use Predefined Method		Show: Files from this instrument	
O Start Favorite			
mth	Name	Remarks Status	
Help Cancel <<< Back	M	AKE YOUR SELECTION	
<b>Obtain Raw Data</b> (Erhålla rådata)	Används för att generera rådata de nödvändiga mätparametrarna	snabbt genom att ställa in a.	
Use Predefined Method (Använd fördefinierad metod)	Används för att utföra mätningar baserat på redan fastställda metoder.		
<b>Start Favorite</b> (Starta Favorit)	Används för att välja en av de va listan med numrerade ikoner.	anligaste metoderna från	

# 6.2 Erhålla rådata

#### 6.2.1 Erhåll rådata med SUNRISE-instrumentet

Välj Obtain Raw Data (Erhålla rådata) och klicka på Next (Nästa) för att öppna dialogrutan Measurement Parameter (Mätparametrar).

I dialogrutan **Measurement Parameter (Mätparametrar)** finns följande alternativ tillgängliga:

Endpoint measurement (Slutpunktsmätningar)	Använd den här knappen för att utföra en enstaka mätning.
Kinetic measurement (Kinetisk mätning)	Denna typ av mätning används för flera mätningar inom ett visst tidsintervall.
<b>Measurement</b> parameters (Mätparameterar)	Klicka på den här knappen för att definiera parametrarna i dialogrutan <b>Measurement Parameter</b> (Mätparametrar)



VARNING DET REKOMMENDERAS STARKT ATT KONTROLLERA ALLA MÄTPARAMETRAR (ÄVEN PARAMETRAR SOM INTE ANVÄNDS I DEN AKTUELLA MÄTNINGEN), EFTERSOM PARAMETRAR FRÅN DEN TIDIGARE MÄTMETODEN FORTFARANDE KOMMER ATT VARA INSTÄLLDA.

#### 6.2.2 Erhåll rådata med INFINITE F50-instrumentet

När ett INFINITE F50-instrument är anslutet visas följande fönster efter att du valt **Obtain Raw Data (Erhålla rådata)**:

tab Were 2	🔷 = Pak	Selection
+ Pate Part of Plate	Plas defoiser (\$R596) - Serier 56 Flat Transport	Absorbance
🕈 Heasurements 🗉	v Patalitation 2	
4 Alantance Alantance 2 Stating 4 Note Raw Concerned to the state of the state		
Hisorilaneous 2	4 × Absoluce	1
iff Connet te lan fingun de lan fingun de lan fingun de lan fingun te lan f	Nacional Editer Series	
Number of plates: 1 (2) Melp Cancel <<< Back	CHOOSE MEASUREMENT P	ARAMETERS

Definiera mätparametrarna enligt listan på vänster sida.

Klicka på **Choose Measurement Parameters (Välj mätparametrar)** För att avsluta definitionen. Se följande instruktioner för **i-control** för ytterligare information.



# 6.3 Use Predefined Method (Använd fördefinierad metod)

Klicka på **Use Predefined Method (Använd fördefinierad metod)** och välj en metod från listan **Filename (Filnamn)**.

De tillgängliga metoderna visas beroende på valet av rullgardinsmenyn **Show** (Visa), dvs. All files (Alla filer), Files from this instrument (Filer från det här instrumentet), My files (Mina filer), Signed files (Signerade filer) eller Last selected methods (Senast valda metoder).

## 6.4 Start Favorite (Starta Favorit)

Klicka på **Start Favorite (Starta Favorit)** för att välja en metod som tidigare definierats som en så kallad favoritmetod som används ofta från listan med numrerade ikoner.

Klicka på **Organize favorites (Organisera favoriter)** för att lägga till eller ta bort metoder från favoritlistan.

#### **Organize Favorites (Organisera favoriter)**

Metodfiler kan dras från listan **Method (Metod)** till en av de 15 ikonerna i listan **Favorite (Favorit)** för att spara den som en favorit.

Organize Fav	orites							
Methods:								
Filename			Remarks					
Mualitative Muanti	Elisa.mth re Elisa.mth ite.mth							Set Remove
Favorite method	ls:							
Empty>	2 <empty></empty>	3 <empty></empty>	4 <empty></empty>	5 <empty></empty>	6 <empty></empty>	< <b>T</b>	<b>8</b> <empty></empty>	
<b>9</b> <empty></empty>	<b>10</b> <empty></empty>	<b>11</b> <empty></empty>	<pre>Empty&gt;</pre>	<b>13</b> <empty></empty>	<b>14</b> <empty></empty>	15 <empty></empty>		
infinite M1000 StartFavorite.	IInstrument Favo mth	orite		Select			ок (	Help Cancel

Dialogrutan **Organize Favorites (Organisera favoriter)** innehåller följande element:

Listan <b>Methods</b> (Metoder)	Listar alla tillgängliga metoder. Filnamnet och motsvarande anmärkningar - om de har angetts - visas.
Knappen <b>Set</b> (Ställ in)	När du har valt en metod från listan <b>Methods Filename</b> (Metoders filnamn) och valt en metodikon klickar du på <b>Set</b> (Ställ in) för att lägga till den valda metoden i <b>Favorites</b> (Favoriter).
Knappen <b>Remove</b> (Ta bort)	Välj en metodikon och klicka på <b>Remove</b> (Ta bort) för att radera en metod från <b>Favorites</b> (Favoriter).

## 6.5 Starta mätning med en fördefinierad eller favoritmetod

Klicka på **Make your selection (Gör ditt val)** respektive **Show Summary (Visa sammanfattning)** i dialogrutan **Select a file (Välj en fil)** för att öppna dialogrutan **Start Measurement (Starta mätning)**. Innan instrumentet börjar mäta måste användaren (om definierat) svara på rapportfrågorna och/eller bekräfta, eller ändra konstanterna.

#### **Dialogrutan Edit Prompts (Redigera prompter)**

Om användarinstruktioner har angetts i metoden visas Edit user prompts (Redigera användarprompt):

Listan Prompt Answer (Promptsvar)	I varje textfält i denna lista kan du skriva in vilken text som helst.
Kryssrutan <b>Req.</b> (Krav)	Kryssrutan <b>Required</b> (Krav) bredvid varje <b>Prompt</b> <b>Answer</b> (Promptsvar)-textfältet fastställer att mätningen endast kan startas när en text har angetts.

Klicka på **Continue (Fortsätt)** för att stänga dialogrutan för prompter. **Fönstret Start Measurement (Starta mätning)** visas. Om den anges i metoden, kommer dialogrutan **Edit Constant (Redigera konstanter)** att visas.

#### Dialogrutan Edit Constants (Redigera konstanter)

Dialogrutan **Edit Constants (Redigera konstanter)** visas endast om den nödvändiga flaggan är inställd för minst en konstant i den valda metoden.

Listan <b>Constants</b> (Konstanter)	I varje textfält i denna lista kan värdet av konstanten eller kommentaren ändras.
Kryssrutan <b>Req.</b> (Krav)	Kryssrutan <b>Req.</b> (Krav) bredvid varje <b>Constant value</b> (Konstant värde)-textfält fastställer att det konstanta värdet måste bekräftas. Gör detta genom att gå vidare till nästa sida.

Om metodkonstanterna har ändrats kan de nya värdena automatiskt överföras och sparas i motsvarande metod. Klicka på **Modify constants in Method (Ändra konstanter i metod)** för att ändra den aktuella metoden med de nya definierade konstanterna. Observera att denna dialogruta endast är tillgänglig för **magellan**användare med användarrättigheterna **Edit methods (Redigera metoder)**.

Klicka på **Continue (Fortsätt)** för att stänga dialogrutan för konstanter. Fönstret **Start Measurement (Starta mätning)** visas.





Anmärkning

Om du använder magellan Tracker eller magellan Standard med användaradministration, kan en standard magellan-operatör (se kapitel 9.6 Användarrättigheter och 9.5.3 Ändra användare) endast redigera de nödvändiga konstanterna. De konstanter som inte har den nödvändiga flaggan visas i grått och kan inte redigeras.

En standard magellan-administratör kan också redigera konstanter utan att den nödvändiga flaggan är inställd.

För en standard magellan programspecialist beror det på vilken metod som används. Med hjälp av en osignerad metod kan användaren redigera konstanter med och utan den nödvändiga flaggan. Med hjälp av en signerad metod kan användaren endast redigera de nödvändiga konstanterna

Efter dessa två valfria sidor visas dialogrutan **Start Measurement (Starta mätning**:

art Measurement			
Measurement			Measurement parameters
Workspace:	06032018-001.wsp	Arb. cycle kin	Measurement mode: Absorbance
Method:	Application Example_en.mth	Modify layout	Measurement wavelength: 492 nm
Sample ID list:		Insert	Read mode: Normal
Instrument			Unit: OD
Use sta	icker	Current: 32.5 °C	
Plate in	Temp. control	Target: 37.0 °C	
Movements.	. Injector control	Optimize Z-position	
Please note:			
	/		
Help Cancel	<<< Back		START
<sup>-</sup> upprutan e <b>asuremen</b> lätning)	t arbetsyta (Automati arbetsyta	Workspace (A enligt definition serad datahan kan bytas ut.	Arbetsyta): Standardfilnamnet på nen i <b>Automated Data Handlin</b> tering) visas. Filnamnet för denr
			<i></i>

Knappen **Arb. cycle kin. ...** (Arbiträr kinetikcykel): Den här knappen är tillgänglig om en kinetisk mätning utförs. Klicka på den här knappen för att visa dialogrutan **Arbitrary Cycle Kinetic** (Arbiträr kinetikcykel). . I den här dialogrutan kan mätningen delas upp i flera delmätningar med olika antal mätcykler och intervalltider. Var och en av dessa delmätningar kan startas när som helst. Arbetsytan kan sparas och mätningen kan återupptas en annan dag.



Anmärkning Detta alternativ är inte tillgängligt för INFINITE F50-instrument.

	Textfältet <b>Method</b> (Metod): Filnamnet för den tidigare valda metoden visas.
	Knappen <b>Modify layout</b> (Ändra layout): Beroende på användarrättigheterna kan plattlayouten ändras. Klicka på knappen <b>Modify layout</b> (Ändra layout) för att öppna dialogrutan <b>Modify layout</b> (Ändra layout). Användaren kan flytta kontroller och ta bort prover, ändra koncentrationen eller ändra mätparametrarna. Ändringarna lagras endast i arbetsytan och inte i den tidigare valda metoden. Den här knappen är inte tillgänglig om du har valt Obtain Raw Data (Erhålla rådata) och efter att du har infogat en prov-ID-lista.
	Textfältet <b>Sample ID List</b> (Prov-ID-lista): Filnamnet för den aktuellt laddade prov-ID-listan visas
	Knappen <b>Insert</b> (Infoga): se även nedan. Klicka på den här knappen för att öppna dialogrutan <b>Insert Sample ID List</b> (Infoga prov-ID-lista). Efter att ha valt prov-ID-lista visas dialogrutan <b>Sample ID List with Plate Layout Preview</b> (Prov-ID-lista med förhandsgranskning av plattlayout) där användaren måste bekräfta valet.
Grupprutan Instrument	Arbetsytorna kommer att lagras med standardnamn för arbetsytorna. När alla plattor har mätts är endast den första arbetsytan öppen. Använd <b>File Open (Öppen fil)</b> eller guiden <b>Evaluate Results</b> (Utvärdera resultat) för att öppna de andra arbetsytorna. <b>Det är inte möjligt att använda</b> <b>staplaren vid kinetiska mätningar.</b>
	Knappen <b>Plate In/Out</b> (Platta in/ut): Klicka för att flytta på platthållaren. Den här knappen ändrar texten beroende på platthållarens position.
	Knappen <b>Movements</b> (Förflyttning): Ytterligare förflyttningar av instrumentet kan hänvisas (filterhållare, spegelhållare osv.)
	Textfältet <b>Current Temperature</b> (Aktuell temperatur): Instrumentets aktuella temperatur visas.
	Textfältet <b>Target Temperature</b> (Måltemperatur): Måltemperaturen, som ställs in i dialogrutan <b>Temperature</b> <b>control</b> (Temperaturkontroll), visas.
	Knappen <b>Temperature Control</b> (Temperaturkontroll): Klicka på den här knappen för att öppna dialogrutan <b>Temperature</b> <b>Control</b> (Temperaturkontroll). Se kapitel 3.1.2 Temperature Control (Temperaturkontroll).
Grupprutan <b>Measurement</b> <b>parameters</b> (Mätparameterar)	I ett textfält listas alla konfigurerade mätparametrar och de kan dubbelkontrolleras.
Grupprutan <b>Please note</b> (Observera)	Viktig information visas. Läs den här informationen och vidta lämpliga åtgärder innan du kör metoden.

En varning visas om layouten definieras för brunnar som inte kommer att mätas.



#### Klicka på Start för att starta mätningen.

Om du kör en metod med flera plattor, visas dialogrutan **Run Measurement (Kör mätning)** efter att du har klickat på **Start**.

lun measurement	
Measurement of plate 1 of 3	
Sample IDs list:	Insert
STOP PAUSE RU	
	Plate in

Den innehåller följande delar:

Measurement of plate (Mätning av platta)	Visar numret på den aktuellt behandlade plattan.
<b>Sample ID List</b>	Infoga lämplig prov-ID-lista för den valda plattan. Prov-ID-
(Prov-ID-lista)	listor måste infogas platta för platta.



Anmärkning
Endast prov-ID-listor som skapats i magellan kan infogas. Import av andra
format stöds inte.

RUN	Klicka på den här knappen för att starta mätningen.
PAUSE	Du når guiden Evaluate Results (Utvärdera resultat) när du klickar på <b>Pause</b> (Paus); ytterligare plattor kan läggas till senare.
STOP	Du når guiden Evaluate Results (Utvärdera resultat) när du klickar på <b>Stop</b> (Stoppa); inga ytterligare plattor kan läggas till.

När mätningen är avslutad visas dialogrutan **Results (Resultat)**, där alla resultat och beräkningar kan visas. Se kapitel 7 Evaluate Results Wizard (Guiden Utvärdera resultat) för mer information.

#### Insert Sample ID List (Infoga en prov-ID-lista)

Alternativet **Insert Sample ID List (Infoga en prov-ID-lista)** används för att ladda in en tidigare definierad prov-ID-lista i en arbetsyta:

Fältet Filename (Filnamn)	Listan visar alla tillgängliga prov-ID-listor (.smp). Välj en prov-ID-lista genom att klicka på ett filnamn.
Fältet <b>Remarks</b> (Anmärkningar)	Tillgängliga kommentarer och anmärkningar visas.



Knappen	Klicka för att öppna standarddialogrutan Windows Open
Import	(Öppna). Externa filer kan importeras. Se kapitel 5.2.3
(Importera)	Importera en prov-ID-lista för mer information.

Välj önskad prov-ID-lista från fältet **Filename (Filnamn)** och klicka på **OK** för att visa prov-ID-listan i fönstret **Sample ID List with Plate Layout Preview (Prov-ID-lista med förhandsgranskning av plattlayout)**. I den här dialogrutan visas listan över prov-ID:n tillsammans med den definierade plattlayouten. Om metoden och prov-ID-listan har olika plattformat visas ett felmeddelande.

När prov-ID-listan väljs i inställningen för den utskrivna rapporten visas den på utskriften.

#### Sample ID List with Plate Layout Preview (Prov-ID-lista med förhandsgranskning av plattlayout) visar prov-ID-listan tillsammans med den definierade plattlayouten. Detta hjälper användaren att se till att rätt prov-ID-lista väljs. Dialogrutan innehåller följande element:

Knappen <b>10%</b>	Klicka för att ändra storleken på den visade plattlayouten så att den visar <b>10 %</b> av plattlayouten på helskärm.
Knappen 100%	Klicka för att ändra storleken på den visade plattlayouten så att den visar <b>100 %</b> av plattlayouten på hela skärmen.
Kryssrutan <b>Zoom</b>	När denna <b>Zoom</b> -funktion är aktiverad kan de nödvändiga brunnarna på plattlayouten markeras och brunnarna förstoras automatiskt så att de passar på hela skärmen.
Kryssrutan Remove unused samples (Ta bort oanvända prover)	Om en endast delvis fylld prov-ID-lista laddas kan användaren radera alla oanvända prov från layouten genom att välja den här kryssrutan. Om en slutlig kontroll (t.ex. i brunn H12) har definierats i layouten flyttas denna kontroll automatiskt till brunnen efter den senast använda provbrunnen. Som standard är alternativet <b>Remove unused samples (Ta bort</b> <b>oanvända prover)</b> vald.

# 6.6 Measurement Status (Mätstatus)

Dialogrutan **Measurement Status (Mätstatus)** visas under instrumentets mätning och innehåller följande element:

Textfältet Status	Textfältet uppdateras under mätningen och innehåller information om mätningens status.
Grupprutan Display in well (Visa i brunn)	Välj hur data ska representeras i brunnarna: Values (Värden): Det uppmätta värdet visas i brunnen. Om mer än ett värde mäts visas alla värden i en brunn.
	<b>Graph</b> (Diagram): Alternativet är aktiverat om ett kinetiskt diagram kan visas. Diagrammet visas så snart mer än ett värde är tillgängligt.
Knappen <b>Pause</b> (Paus)	Om du vill pausa en kinetisk mätning vid nästa möjliga tidpunkt klickar du på knappen <b>Pause</b> (Paus). I dialogrutan <b>Pausing Kinetic</b> (Pausa kinetik) klickar du på knappen <b>Plate out</b> / <b>Plate in</b> (Platta ut/in) för att flytta platthållaren ut ur eller in i instrumentet. Klicka på <b>Continue</b> (Fortsätt) för att fortsätta med den kinetiska omgången.
Knappen <b>Break (Bryt)</b>	Klicka på den här knappen för att avbryta mätningen vid nästa möjliga tidpunkt.



#### 6. Start Measurement Wizard (Guiden starta mätning)

Displayen <b>Plate View</b> (Plattvy)	Displayen <b>Plate View</b> (Plattvy) visar en schematisk bild av mikroplattan. Enligt alternativet <b>Display in wells</b> (Visa i brunnar) anpassas innehållet i brunnarna. Brunnarna uppdateras så snabbt som instrumentet levererar data. När injektion eller dispensering utförs markeras den aktuellt använda brunnen automatiskt med en annan färg.
Textfältet <b>Cycle</b> (Cykel)	Om en kinetisk mätning utförs, visar textfältet <b>Cycle</b> (Cykel) det aktuella cykelnumret och det maximala antalet kinetikcykler.
Displayen <b>Graph</b> (Diagram)	Om en kinetisk mätning eller en mätning med multipla inmärkningar utförs, visas det kinetiska diagrammet eller diagrammet med multipla inmärkningar för den aktuellt valda brunnen i en större storlek på höger sida av dialogrutan <b>Measurement Status</b> (Mätstatus).
Grupprutan <b>Previous cycle</b> (Föregående cykel)	Så länge överföringen av data för den aktuella cykeln inte påbörjades innehåller textfälten <b>Raw data</b> (Rådata) och <b>Temperature</b> (Temperatur) värdet samt temperaturen för den föregående uppmätta cykeln av den aktuellt valda brunnen.
Grupprutan <b>Time</b> (Tid)	Textfältet <b>Elapsed time</b> (Förfluten tid) visar den tid som förflutit sedan mätningen inleddes.
	Textfältet <b>Exp. run time</b> (Förväntad körtid) visar den förväntade tidsperioden för hela mätningen.

Skalning av y-axeln i diagrammet är möjlig genom att välja antingen Auto select range (Välj intervall automatiskt) eller Select range (Välj intervall) (MIN/MAX).



Anmärkning Det är möjligt att välja flera brunnar på plattan för att visa de valda kinetiska kurvorna i ett diagram. Vid kinetiska mätningar sparas de hämtade data automatiskt var 30:e minut i magellans loggfilskatalog (se

kapitel 3.3 Log Files (Loggfiler)).

2023-07

# 7. Evaluate Results Wizard (Guiden Utvärdera resultat)

# 7.1 Inledning

Guiden **Evaluate Results (Utvärdera resultat)** är utformad för att hjälpa användaren att analysera mätresultat. Rådata, utvärderingsdata och utvärderingsparametrar kan visas och data kan utvärderas på nytt.

#### Sammanfattning av arbetsflöde

Starta guiden **Evaluate Results (Utvärdera resultat)** och välj en befintlig arbetsyta för utvärdering (Dialogrutan **Select a File (Välj en fil)**)

Klicka på **Make your selection (Gör ditt val)** för att öppna dialogrutan **Results** (**Resultat**) där resultaten kan visas eller skrivs ut (Fliken **Evaluate Results** (**Utvärdera resultat**)), och vid behov kan metoden som används ändras (Fliken **Edit Method (Redigera metod)**).

Klicka på **Finish (Slutför)** och använd dialogrutan **Save (Spara)** för att spara en arbetsyta som innehåller metoddefinitioner, instrumentdata och en prov-ID-lista om konfigurerad.

Utvärderade data lagras inte i arbetsytan utan beräknas på nytt varje gång arbetsytan öppnas.

# 7.2 Select a File (Välj en fil)

Dialogrutan Select a File (Välj en fil) innehåller följande element:

Listan <b>Filename</b> (Filnamn)	Listan <b>Filename</b> (Filnamn) innehåller en lista över alla filer (och deras kommentarer) i standardarbetsytans katalog.
Kombinationsrutan <b>Show</b> (Visa)	I Kombinationsrutan <b>Show</b> (Visa) kan den visade listan med filer ändras enligt valet:
	<ul> <li>Files from this instrument (Filer från det här instrumentet)</li> </ul>
	<ul> <li>My files (Mina filer): Det här alternativet är tillgängligt om användaradministrationen är aktiverad (den är alltid aktiverad i magellan Tracker).</li> </ul>
	<ul> <li>Signed files (Signerade filer): endast tillgängliga för magellan Tracker</li> </ul>
	<ul> <li>Anpassade definitioner är möjliga genom att klicka på knappen; se File Selection Criteria (Kriterier för filval) nedan.</li> </ul>
	• Example files (Exempelfiler): endast tillgängliga om de har installerats.



#### 7.2.1 Kriterier för filval

File Selection Criteria	
Workspace criteria <ul> <li>All files</li> <li>Files from this instrument</li> <li>My files</li> <li>Signed files</li> <li>Containing</li> </ul>	OK Cancel
Workspace file created/modified No time limitation Today Last week	
O Last D days	Help

Klicka på knappen för att öppna dialogrutan File Selection Criteria (Kriterier för filval).

Välj ett av alternativen under Workspace Criteria (Kriterier för arbetsyta) eller ange tecken i textfältet Containing (Innehåller) som måste finnas i Filename (Filnamn). Ett tidsfilter kan tillämpas genom att välja ett av alternativen under Workspace file created/modified (Arbetsyta skapad/ändrad) eller ange ett nummer i fältet Last ... days (Senaste ... dagar). Klicka på OK för att spara inställningarna.

Välj den arbetsyta som ska utvärderas i listan Filename (Filnamn) och klicka på Make your selection (Gör ditt val).



#### Anmärkning

Om ett SUNRISE-instrument är anslutet, men en INFINITE F50-metod är vald, visas dialogrutan Instrument Mismatch (Missmatchade instrument):

• Anslut till rätt instrument.



# 7.3 Fönstret Workspace Overview (Översikt över arbetsyta)

Fönstret **Workspace Overview (Översikt över arbetsyta)** visar alla tillgängliga data för den valda arbetsytan.

Om den användare som arbetar för tillfället har rätt **magellan**-rättigheter, är det också möjligt att göra vissa ändringar i metoden genom att klicka på fliken **Edit Method (Redigera metod)** och göra en ny beräkning av de uppmätta data.



Workspace Overview (Översikt över arbetsytan) består av följande element:

- Fönstret Plate layout (Plattlayout): utvalda data för varje brunn visas i plattlayouten. Om en arbetsyta med flera plattor öppnas visas en "räknare" för att växla mellan de olika plattorna i det övre högra hörnet.
- Verktygsfält: för att redigera, zooma, växla mellan kinetikcykler och att visa sammanfattningen av en vald brunn.
- Fönstret **Data Information (Datainformation)**: data som inte kan tilldelas en enstaka brunn visas i textrutan (t.ex. mätparametrar, definition av cutoff, resultatstatistik osv.).
- Kontrollfältet i fliken Evaluate Results (Utvärdera resultat): alla tillgängliga namn på dataset visas. Om du väljer ett dataset visas antingen data i plattlayouten eller data i fönstret Data Information (Datainformation).
- Kontrollfältet i fliken Edit Method (Redigera metod): inställningarna och parametrarna för den laddade metoden visas.

Storleken på de tre sistnämnda fönstren kan ändras med hjälp av de delningslister som skiljer fönstren åt.



#### Plattlayout

Vyn visar en schematisk layout av en mikroplatta. Raderna markeras i alfabetisk ordning, kolumnerna i numerisk ordning.

I kontrollfältet på vänster sida visas alla tillgängliga data. De utvalda data visas i plattans layout, i textpanelen eller i ett separat diagramfönster, beroende på vilken typ av utvalda data.

#### Sammanhangsberoende meny för plattlayouten

Om du vill visa en sammanfattning av alla data för en brunn, markera brunnen och välj **Summary (Sammanfattning)** från menyn med höger musknapp.

#### Knappar i verktygsfältet

De tillgängliga knapparna i verktygsfältet beskrivs i följande punkter.

Beroende på vilka data som valts visas olika knappar:

när Instrument data (Instrumentdata) väljs, bli knappen Edit (Redigera) tillgänglig;

när Concentration (Koncentration) eller Kinetiska data väljs, blir knappen Curve/Graph (Kurva/diagram) tillgänglig.

Följande knappar är endast tillgängliga i guiden **Evaluate results (Utvärdera resultat)**.

Edit (Redigera)	Klicka på <b>Edit (Redigera)</b> för att redigera eller dölja rådata.
Knappen 🛄	Klicka på den här knappen för att visa dialogrutan <b>Summary (Sammanfattning)</b> .
Knappen 10%	Klicka på den här knappen för att ändra storleken på den visade plattlayouten så att den visar <b>10 %</b> av plattlayouten på helskärm.
Knappen 100%	Klicka på den här knappen för att ändra storleken på den visade plattlayouten så att den visar <b>100 %</b> av plattlayouten på helskärmen.
Kryssrutan Zoom	När kryssrutan <b>Zoom</b> är markerad kan de nödvändiga brunnarna på plattlayouten markeras och de märkta brunnarna förstoras automatiskt så att de passar hela skärmen.
Rullgardinsmenyn Measurement cycle (Mätcykel)	Om mätningen är en kinetisk mätning kan önskad mätcykel väljas i rullgardinsmenyn.

#### Fönstret Data Information (Datainformation)

Du kan till exempel klicka på **Miscellaneous (Övriga)** data i fliken **Evaluate Results (Utvärdera resultat)** för att visa motsvarande information i fönstret **Data Information (Datainformation)**.



# 7.4 Fliken Evaluate Results (Utvärdera resultat)

Alla tillgängliga namn på dataset visas i kontrollfältet **Available Data** (Tillgängliga data). Det finns flera olika visningsalternativ:

- Utvalda data som visar ett värde för varje brunn visas i fönstret för plattlayout. Den kan visas som värden eller som pseudofärger tillsammans med en färgskala.
- Kinetiska data kan visas som ett litet diagram i varje brunn.
- Utvalda data som utgör en datainsamling och som inte kan tilldelas en enstaka brunn visas i fönstret för datainformation (t.ex. QC-valideringskriterier).
- Om ett diagram (standardkurva, kinetik, enzymkinetik, spädningsserier, multipla inmärkningar, spektra) väljs, visas dialogrutan **Graph (Diagram)**.

Om en mätsekvens som består av individuella oberoende mätningar har utförts visas resultaten av varje mätning när det relevanta mätningsnumret väljs från en rullgardinsmeny i verktygsfältet (t.ex. kinetikcykler i en kinetisk omgång).

Vid behov kan mätresultaten ändras genom att klicka på knappen **Edit** (**Redigera**). Manipulerade data är markerade med ett ~-tecken.



#### VARNING

#### DENNA FUNKTION ÄR ENDAST AVSEDD FÖR FORSKNINGSÄNDAMÅL OCH FÅR ALDRIG ANVÄNDAS FÖR DIAGNOSTISKA TESTER.

I varje brunn av plattlayouten kan tre rader av data visas. Beroende på vilka kriterier som valts i kontrollpanelen i det vänstra fönstret, visas olika värden.

#### 7.4.1 Menyer

<b>File</b> (Fil)	ASCII File Export (Export av ASCII-fil) ASTM File Export (LIS) (Export av ASTM - LIS) Sample Tracking Export (Spårning av prov) Excel Export (Export av Excel) Method Export (Export av metod) Print (Skriv ut) Print Preview (Förhandsgranskning av utskrift) Printer Setup (Skrivarinställningar) Printout Font (Teckensnitt för utskrift) Plate to Plate QC (QC platta-till-platta)
<b>Edit</b> (Redigera)	Copy to Excel (Kopiera till Excel) Copy to ASCII Format (Kopiera till ASCII-format) Paste from ASCII Format (Klistra in från ASCII-format) Insert Sample ID List (Infoga en prov-ID-lista) Recalculate with another Method (Beräkna på nytt med en annan metod)
Instrument	Movements… (Förflyttningar) Temperature Control… (Temperaturkontroll) Start Measurement… (Starta mätning)

## 7.4.2 Verktygsfältet: File (Fil)

#### ASCII File Export (Export av ASCII-fil)

Klicka på **ASCII File Export (Export av ASCII-fil)** för att exportera data som ASCII-filer (.asc).

Innan exportproceduren inleds måste data väljas i dialogrutan **Data Export** (Dataexport) via fliken Edit method  $\rightarrow$  Data handling  $\rightarrow$  Data export.

Välj den mapp där exporterade data ska lagras. Ange ett filnamn (.asc).

När du klickar på **Save (Spara)** i dialogrutan **Save as (Spara som)** kommer de dataset som anges i dialogrutan **Data export (Dataexport)** att sparas som en ASCII-fil.

#### ASTM File Export (LIS) (Export av ASCTM-fil - LIS)

I menyn **File (Fil)**, klicka på **ASTM File Export (Export av ASTM-fil)** för att exportera data i ASTM-format till kompatibla laboratorieinformationssystem.

Välj de data som ska exporteras från dialogrutan **Data Export (Dataexport)** via fliken **Edit method**  $\rightarrow$  **Data handling**  $\rightarrow$  **Data export**.

#### Sample Tracking Export (Export av spårning av prov)

I menyn File (Fil), klicka på Sample Tracking Export (Export av spårning av prov) för att exportera data till Sample Tracking-systemet (Spårning av prov).

Välj de data som ska exporteras från dialogrutan **Data Export (Dataexport)** via fliken **Edit method**  $\rightarrow$  **Data handling**  $\rightarrow$  **Data export**.

#### **Excel Export (Export av Excel)**

Klicka på Excel Export (Export av Excel) för att exportera data till Excel.

Välj de data som ska exporteras från dialogrutan **Data Export (Dataexport)** via fliken **Edit method**  $\rightarrow$  **Data handling**  $\rightarrow$  **Data export**.

#### Metod Export (Export av metod)

#### Klicka på Method Export (Export av metod) för att exportera en metod.

Välj den mapp där metoden ska sparas och ange önskat filnamn. Dessutom kan några kommentarer läggas till i textfältet **Remarks (Anmärkningar)**.

#### Print (Skriv ut)

I menyn **File (Fil)**, klicka på **Print... (Skriv ut)** eller tryck på **SHIFT-P** för att skriva ut data enligt definitionen i inställningen **Printed Report (Utskriven rapport)** från metoden. Standarddialogrutan Windows **Print... (Skriv ut)** öppnas. Där kan du välja skrivare, sidintervall och antal kopior.

#### Print Preview (Förhandsgranskning av utskrift)

Klicka på **Print Preview... (Förhandsgranskning av utskrift)** för att se exakt hur en utskrift kommer att se ut utan att behöva skriva ut dokumentet.

#### Printer Setup (Skrivarinställningar)

Klicka på **Printer Setup... (Skrivarinställningar)** för att ändra den aktuella skrivaren eller ändra skrivarinställningarna.

Standarddialogrutan Windows **Print Setup (Utskriftsinställning)** visas. Välj skrivare, pappersstorlek och riktning för utskriften.



#### Printout Font (Teckensnitt för utskrift)

Klicka på **Printout Font... (Teckensnitt för utskrift)** för att välja ett visst teckensnitt, stil, storlek, färg, osv. för framtida utskrifter.

Dessa inställningar har ingen effekt på skärmvisningen. Stora teckensnitt bör undvikas för att undvika fel eller felaktig information av data.

#### QC platta-till-platta

När **Plate to Plate QC (QC platta-till-platta)** konfigureras i metoddefinitionen, jämför **magellan** det aktuellt observerade medelvärdet för en kontroll mot det medelvärde och den standardavvikelse som tidigare definierats för den här metoden. (Se kapitel 4 Create/Edit a Method Wizard (Guiden Skapa/redigera en metod) ).

#### Levy-Jennings-diagram

Klicka på Plate to Plate QC (QC platta-till-platta) och Levy-Jenningsdiagrammet visas. Levy-Jennings-diagrammet visar trenden för kontrollvärdena.



På y-axlarna tillämpas de datavärden (rådata, koncentrationer osv.) som anges i metoddefinitionen på tidsskalan dagar (x-axlar).

Kontrollvärdena visas tillsammans med medelvärdet och standardavvikelsernas avståndslinjer (1s, 2s, 3s).

När markören placeras på en mätpunkt visas ytterligare information i verktygstips (arbetsytans namn, osv.). Använd kryssrutorna för de enstaka elementen i listan **Show (Visa)** för att välja olika kontroller som ska visas.

#### Time Range (Tidsintervall)

Klicka på **Time Range (Tidsintervall)** för att utvärdera arbetsytan inom ett definierat tidsintervall.

#### Select Workspaces (Välj arbetsytor)

Klicka på **En-/Disable (Avaktivera/inaktivera)** för att aktivera eller inaktivera vissa arbetsytor för utvärdering.



#### 7. Evaluate Results Wizard (Guiden Utvärdera resultat)

Enable/Disable workspaces
<ul> <li>✓ 20070521-001.wsp</li> <li>✓ 20070521-002.wsp</li> <li>✓ 20070521-003.wsp</li> <li>✓ 20070521-004.wsp</li> <li>✓ 20070521-005.wsp</li> <li>✓ 20070521-006.wsp</li> </ul>
OK Cancel Help

#### Datablad

Klicka på View summary (Visa sammanfattning) för att öppna dialogrutan Data sheet (Datablad):

Kolumnen <b>Workspace</b> (Arbetsyta)	En lista över de utvärderade arbetsytorna. Till vänster om arbetsytans namn visar ett grönt kryss uppfyllelse, ett rött kryss på misslyckande i utvärderingen av Westgard <sup>®</sup> -reglerna (se användarhandboken för <b>magellan</b> för mer information om <b>Westgard-regler</b> ).
Kolumnen Control (Kontroll)	De olika kontrollerna listas.
Kolumnen <b>Value (Värde)</b>	Medelvärdet för varje kontroll visas.
Kolumnerna Westgard® rule evaluation (Utvärdering av Westgard-regel)	De individuella Westgard®-reglerna utvärderas med <b>OK</b> för en kontrollregel som inte har brutits eller <b>x</b> för en kontrollregel som har brutits enligt tabellen nedan (se Westgard-regler nedan).
Fältet Comments (Kommentarer)	Det är möjligt att skriva kommentarer om utvärderingen.
Knappen Print Preview (Förhandsgranskning av utskrift)	Förhandsgranskning av utskrift med Levy-Jennings- diagrammet, kontroll av Westgard-regler och kommentarer.



#### Westgard-regler

Westgard®-reglerna är en QC-modell med flera regler:



#### 7.4.3 Verktygsfältet: Edit (Redigera)

Menyn Edit (Redigera) innehåller följande val:

#### Copy to Excel (Kopiera till Excel)

Kommandot **Copy to Excel (Kopiera till Excel)** tillåter användaren att exportera data direkt till ett MS Excel-kalkylblad.

#### Copy in ASCII-Format (Kopiera i ASCII-format)

Välj **Copy in ASCII-Format (Kopiera i ASCII-format)** för att kopiera de data som finns i de valda brunnarna i plattlayouten till urklippet.

Denna funktion gör det möjligt att överföra brunnsdata till andra programvaror.

#### Paste from ASCII-Format (Klistra in från ASCII-format)

I menyn Edit (Redigera) klickar du på Paste (Klistra in) eller tryck på CTRL-V.

Innehållet i utvalda data kommer att klistras in från urklippet med ASCIIdataformatet. Funktionen gör det möjligt att överföra data från andra programvaror (t.ex: Excel) i brunnarna.

Först måste de relevanta data väljas med musen i den andra programvaran. Därefter måste de utvalda data kopieras och överföras till urklippet. Raderna måste skiljas åt med radmatning och kolumnerna med tabstoppar. Detta görs automatiskt i Excel när du markerar flera celler.

Funktionen **Paste (Klistra in)** klistrar in de kopierade data i ASCII-format från urklippet i brunnarna. Data infogas alltid från position A1. Om data inte är numeriska sätts värdet till 0.

Vid händelse av oavslutad kinetik med definierat interval kommer en fråga att visas. Här bestämmer du om de utvalda data ska skriva över de ursprungliga data eller om data ska bifogas. Tidsintervallet hämtas från mätparametrarna.

#### Insert Sample ID List (Infoga en prov-ID-lista)

Detaljerad information finns på 6.5 Starta mätning - Insert Sample ID List (Infoga en prov-ID-lista).

#### Recalculate with another Method (Beräkna på nytt med en annan metod)

En ny beräkning baserad på inställningarna för en nyvald metod kommer att utföras.

Om ett fel uppstår under den nya beräkningen stoppas proceduren.

Klicka på det här alternativet för att öppna dialogrutan **File open (Öppen fil)**. I en ytterligare **Remarks (Anmärkningar)**-fält visas en textbeskrivning av filen - om den har angetts (endast tillgänglig i **magellan** Tracker). Metoden måste väljas antingen genom att dubbelklicka på metodfilen eller genom att markera metodfilen och klicka på knappen **Open (Öppna)**.

När du har valt metod, visas dialogrutan **Calculating... (Beräkna)**. Den här dialogrutan är endast till för visning och innehåller inga element för redigering. Den stängs när beräkningen är klar.





#### 7.4.4 Verktygsfältet: Instrument

För detaljerad information om **Movements (Förflyttningar)** och **Temperature Control (Temperaturkontroll)** se kapitel 3.1 Instrumentalternativ.

#### Starta mätning

Med det här alternativet kan mätningen startas på nytt med den aktuellt laddade metoden och aktuella data skrivs över om **YES (Ja)** har valts i dialogrutan **magellan**.

Om den kinetiska omgången är ofullständig kan cykler läggas till (inte tillgängligt för instrument i Infinite-serien).

För detaljerad information se 6.5 Starta mätning med en fördefinierad eller favoritmetod.

#### 7.4.5 Fönstret plattlayout

De data som ska visas i en brunn när du öppnar en arbetsyta kan ställas in när du definierar metoden (Automated data handling  $\rightarrow$  view results after measurement  $\rightarrow$  More...).

I varje enstaka brunn syns tre rader, och följande data visas som standard:

1:a raden: layout

2:a raden: replikatinformation

3:e raden: På den tredje raden används de konventioner som beskrivs i följande kapitel, beroende på vilka data som valts (se 7.4.6 Särskilda karaktärer).



#### 7.4.6 Särskilda karaktärer

"None" (Ingen)	Inga data visas.
"~" tilde	En tilde läggs till för det manipulerade värdet av en brunn (simulerad, redigerad), till exempel: 0.354~
"( )" parentes	Parenteser används för det dolda värdet för en brunn, till exempel: (0.354)
"!" utropstecken	Ett utropstecken läggs till för det eliminerade värdet av en brunn; till exempel: 1.0.354
"#" fyrkant	Koncentrationsvärdena markeras med ett "#" när den beräknade koncentrationen ligger utanför standardkurvans intervall, vilket innebär att koncentrationsvärdet har extrapolerats. Till exempel: #13.75
" * " asterisk	En asterisk markerar värden som har uppmätts med alternativet "Använd förstärkningsreglering", som korrigerar (= sänker) förstärkningen.
Motsvarande felmeddelande visas.	Medelvärdet för en brunn är oanvändbart (t.ex.: "lamp low", division med noll, negativ logaritm).
Ingen färg, den tredje raden är tom	Ingen formel tillämpas på brunnen eller Replikatbrunn om "average s" (medelvärde s) eller "v calculation" (v beräkning) eller Inga cutoff-resultat.
Data från den första replikatet döljs eller elimineras	Genomsnittligt värdet erhålls utan att använda dolda data, men det genomsnittliga värdet visas i brunnen för det första replikatet
"< MIN"	Den beräknade koncentrationen ligger under den lägsta standarden.
"MltPt"	Flera punkter; en unik koncentration kan inte beräknas.
"> MAX"	Den beräknade koncentrationen ligger över den högsta standarden.
"NoCalc"	Indata har inte hittats eller beräkningsfel; beräkningen är inte möjlig

#### Genvägar

När du visar transformerade data kan du trycka på CTRL-SHIFT för att visa den definierade formeln på den andra raden så länge du trycker på tangenten.

#### 7.4.7 Kontrollfält Instrumentdata

Användaren kan välja rådata, spektrumdata osv., och deras statistik i enlighet med den utförda mätningen.

#### 7.4.8 Kontrollfält Reducerade data

Användaren kan välja reducerade data, t.ex. differensdata enligt den utförda mätningen.



#### 7.4.9 Kontrollfält Transformerad data

Användaren kan välja den transformation som krävs för att visa de beräknade resultaten, inklusive statistik, färgad display osv.

#### 7.4.10 Kontrollfält Kinetiska parametrar

Användaren kan visa de kinetiska utvärderingsparametrarna, t.ex. medel/max lutning, starttid osv., enligt mätdefinitionen.



Anmärkning Oanvändbara data (t.ex. överflödsvärden) ignoreras vid beräkning av kinetiska data.

#### Definiera reducering av kinetiska data i brunn

Välj **Edit kinetic settings... (Redigera kinetiska inställningar)** från den högra musens sammanhangsberoende meny när en brunn är vald. Dialogrutan för reducering av kinetiska data visas (se kapitel 4.3.6 Kinetik: Reducering av kinetiska data för mer information).

För den valda brunnen kan särskilda parametrar definieras. De kan också kopieras till andra brunnar (välj **Copy kinetic settings (Kopiera kinetikinställningar)** från högra musens snabbmeny, klicka på brunnen/brunnarna där inställningarna ska kopieras och välj **Paste kinetic settings... (Klistra in kinetiska inställningar)**).

Om du ändrar inställningarna för reducering av kinetisk data på fliken Method (Metod) kommer de brunnsspecifika definitionerna att skrivas över.

#### Dialogrutan Graph (Diagram): Enzyme kinetics (Enzymkinetik)

Dialogrutan **Enzyme Kinetics Graph (Diagram för enzymkinetik)** tillåter användaren att visa diagrammet för enzymkinetik.

# Sammanhangsberoende meny av Enzyme Kinetics Graph (Diagram för enzymkinetik)

Genom att högerklicka på diagrammet visas en sammanhangsberoende meny.

Hårkorsmarkör	En hårkorsmarkör visas för att hjälpa till att placera markören på särskilda punkter i diagrammet.
Kopiera till Urklipp	Kopierar diagrammet till urklippen som en bitmapp; bitmappen kan överföras till alla Windows-program med hjälp av funktionen klistra in.
Save as bitmap (Spara som bitmapp)	Sparar diagram som en bitmappfil.
Print (Skriv ut)	Skriver ut diagrammet.
Print… (Skriv ut) Printer Setup… (Skrivarinställningar)	Skriver ut diagrammet. Visar dialogrutan <b>Printer Setup</b> (Skrivarinställningar) där skrivarinställningarna kan definieras.



**TECAN**.

#### 7.4.11 Kontrollfält Koncentrationer

Välj enstaka koncentration, medel eller genomsnittlig enstaka koncentration för att visa resultaten, beräknade enligt standardkurvan:

- Single conc. (Unit)
   Koncentrationen för ett individuellt replikat.
- Mean conc. (Unit) Medelvärdet av replikat används för beräkning av koncentrationen. Om replikat med olika spädningar definieras kommer medelkoncentrationen inte att vara tillgänglig.
- Average single conc. (Unit) Koncentrationen beräknas för varje replikat. Därefter beräknas medelvärdet av koncentrationerna.
- Ytterligare koncentrationer (enstaka koncentration, medelkoncentration och genomsnittlig koncentration) om de finns tillgängliga.
- Diagram: Standard Curve (Standardkurva) Klicka på det här alternativet för att öppna dialogrutan Graph: Standard Curve (Diagram: Standardkurva)
- Intercepts (Skärningspunkter):
- Värden för koncentrationstransformationer, om de definieras
- IC50, r-IC50, Diagram: spädningsserier

#### Dialogrutan Graph: Standard Curve (Diagram: Standardkurva)

Öppna den här dialogrutan via kontrollfältet eller genom att klicka på Curve (Kurva) i verktygsfältet för att visa och redigera Standard Curve (Standardkurva).

Den här dialogrutan har en egen meny med alternativ för att spara standardkurvan, ändra analystyp eller jämföra ett antal kurvor. Diagram kan exporteras som en bitmapp som sedan kan inkorporeras i dokument i andra programvaror.

# **Dialogrutan Graph: Standard Curve (Diagram: Standardkurva)** innehåller följande element:

#### Filmeny

Filmenyn innehåller följande kommandon:

 New (Ny) väljs för att skapa en ny standardkurva, X- och Y-värdena kan definieras i en dialogruta.
 Ange uppmätta värden (Y-värde) och koncentration (X-värde). Om alternativet Exclude (Exkludera) väljs, kommer motsvarande punkt inte att tas med i beräkningen av diagrammet.

När du är klar klickar du på knappen OK.



- **Open** (Öppna) väljs för att öppna en tidigare sparad standardkurva. En standardkurva som sparats i filformatet .std kan öppnas. När detta kommando utförs visas diagrammet tillsammans med de redan visade standardkurvorna.
- Close (Stäng) Om flera kurvor är öppna kan individuella kurvor stängas med detta alternativ.

Den aktuella kurvan måste väljas från den tillgängliga listan och stängas genom att klicka på knappen **OK**.

- Save/Save as (Spara/Spara som) väljs för att spara en standardkurva för ytterligare utvärderingar.
   En standardkurva kan sparas som en .std fil. Filen kan inkluderas i metoden genom att växla till fliken Data. Om användaren har valt Save (Spara) och kurvan tidigare har sparats, sparas den utan att några prompter visas. Om det inte är fallet uppmanas användaren att ange ett namn för den nya kurvan.
- **Export** (Exportera) väljs för att exportera en standardkurva som en bitmappfil.
- **Print...** (Skriv ut) väljs för att skriva ut standardkurvan. Detta kommando skriver ut det aktuella diagrammet.
- Printer Setup... (Skrivarinställningar)
   Om du väljer det här alternativet visas dialogrutan Printer Setup... (Skrivarinställningar).
- Exit (Avsluta) väljs för att avsluta dialogrutan för standardkurvan. Om några ändringar har gjorts, t.ex. en ändring av interpoleringsprocessen, görs en ny beräkning.



Anmärkning Standardkurvor som läggs till med New (Ny) eller Open (Öppen) är endast till för jämförelse. När du stänger dialogrutan Standardkurva tas kurvorna bort.

# **EXPERTENS KUNSKAP**

En sparad fil med standardkurva kan användas i metoder som inte har standarder i layouten för att beräkna provkoncentrationer. Se 4.3.8 Koncentrationer:Standardkurva, **Standarder från extern fil**.

#### Redigeringsmenyn

Menyn Edit (Redigera) innehåller följande kommandon:

- Välj Copy (Kopiera) för att kopiera diagrammet till urklippet som en bitmapp, som kan överföras till alla Windows-program med hjälp av funktionen kopiera/klistra in.
- Välj Properties (Egenskaper) för att ändra analystyp, axlar, titel osv.

• Välj **Standard Curve** (Standardkurva) för att öppna dialogrutan Standardkurva. I den här dialogrutan kan standardpunkter uteslutas eller inkluderas.

När du väljer den relevanta kurvan i den tillgängliga listan och klickar på **OK** kommer alla baspunkter att visas i fönstret som visas. Aktivera eller uteslut en punkt genom att klicka på alternativet **Exclude** (Exkludera). Endast de punkter som inte har markerats som uteslutna kommer att användas vid beräkningen av standardkurvan.

Endast baspunkter som har erhållits från rådata, som har matats in manuellt eller som har laddats in ytterligare kan ändras. Baspunkter som erhålls från transformation, till exempel, kan bara uteslutas.

 Conc. Range (Koncentrationsintervall): Dialogrutan Concentration Range (Koncentrationsintervall) visas. Välj mellan Display all (Visa alla) och Display range (Visa intervall). Om Display range (Visa intervall) har valts och Min.- och Max.-gränser definieras, visas endast koncentrationsvärden inom det angivna intervallet i fönstret plattlayout.

#### Menyn View (Visa)

Menyn View (Visa) innehåller följande kommandon:

- Välj **Audit trail** för att visa audit trail för en extern standardkurva. Detta alternativ är endast tillgängligt för **magellan** Tracker.
- Välj Statistics (Statistik) för att visa statistiken över de erhållna värdena.
   För att bedöma om någon drift eller förändring har skett med läsaren under en tidsperiod är det möjligt att jämföra en serie standardkurvor med hjälp av statistiska värden.

Om ett antal kurvor laddas beräknas medelvärdet, standardavvikelsen och variationskoefficienten för var och ett av dem.

- Intercepts... (Skärningspunkter) öppnar dialogrutan Intercepts (Skärningspunkter). Dialogrutan Intercepts (Skärningspunkter) visar resultatet av de definierade värdena för skärningspunkter.
- Average Standard Curve (Genomsnittlig standardkurva) visar standardkurvan som ett medelvärde för de utförda försöksgrupperna. Det här alternativet kan bara aktiveras om det finns ett antal kurvor tillgängliga i fönstret. En genomsnittlig kurva genereras och visas med hjälp av insamlade data av kurvor.

#### Menyn Help (Hjälp)

Välj menyn Help (Hjälp) för att öppna dialogrutan Help (Hjälp) i magellan.

Click on points (Klicka på punkter)

**Click on points (Klicka på punkter)** kommer att dölja/visa dem i standardkurvan. När en punkt har dolts justeras linjen i diagrammet automatiskt i enlighet med detta och punkten representeras som en genomskinlig symbol i stället för en solid symbol.

Tips: Detta kan ångras genom att trycka på CTRL+Z.

#### X- och Y-markörens verktygstips

När markören inte flyttas under en kort stund, visas ett verktygstips med X- och Y-koordinaterna för markörens aktuella position.



# **EXPERTENS KUNSKAP**

Om fler än **sex** standardkurvor visas i diagrammet, visar endast en liten förklaring på den högra sidan, diagrammets etikett. Om du vill visa anpassningsfunktionerna respektive parametrarna för kurvanpassning, välj



Standard Curve Data (Data för standardkurva) från den sammanhangsberoende menyn (se nedan).

#### Sammanhangsberoende meny för ett diagram för standardkurva

Genom att högerklicka på diagrammet visas en sammanhangsberoende meny.

Hårkorsmarkör	En hårkorsmarkör visas för att hjälpa till att placera markören på särskilda punkter i diagrammet.
Single points (Enstaka punkter)	Visa replikatvärden i stället för medelvärden.
Copy to Clipboard (Kopiera till Urklipp)	Kopierar diagrammet till urklippet som en bitmapp, som kan överföras till alla Windows-program med hjälp av funktionen klistra in.
Save as Bitmap (Spara som bitmapp)	Sparar diagrammet som en bitmappfil.
Print… (Skriv ut)	Skriver ut diagrammet.
Printer Setup (Skrivarinställning ar)	Visar dialogrutan <b>Printer Setup</b> (Skrivarinställningar).
Undo changes (Ångra ändringar)	Avbryter nyligen genomförda åtgärder.
Zoom to 100% (Zooma till 100 %)	Ställer tillbaka diagramvisningen till 100 % om zoomfaktorn har ändrats genom att välja rektangulära områden i diagrammet.
Standard Curve Data (Data för standardkurva)	Visar parametrarna för kurvanpassning för den/de visade standardkurvan/standardkurvorna.
Intercepts (Skärnings- punkter)	Öppnar dialogrutan <b>Intercepts (Skärningspunkter)</b> som visar resultatet av de definierade värdena för skärningspunkterna.
Properties (Egenskaper)	Ändrar egenskaperna för diagramvisningen. Diagrammets titel, axelformat, kurvstil osv., kan anpassas.
Help (Hjälp)	Öppnar standarddialogrutan hjälp i magellan.

#### Dialogrutan Graph: Dilution Series (Diagram: Spädningsserie)

Dialogrutan Graph: Dilution Series (Diagram: Spädningsserie) visar spädningsdiagrammet med den angivna skärningspunkten.

#### Sammanhangsberoende meny för spädningsdiagram

Genom att högerklicka på diagrammet visas en sammanhangsberoende meny.

Hårkorsmarkör	En hårkorsmarkör visas för att hjälpa till att placera markören på särskilda punkter i diagrammet.
Copy to Clipboard (Kopiera till Urklipp)	Kopierar diagrammet till urklippet som en bitmapp; bitmappen kan överföras till alla Windows-program med hjälp av funktionen klistra in.
Save as bitmap (Spara som bitmapp)	Sparar diagrammet som en bitmappfil.

Print (Skriv ut)	Öppnar dialogrutan <b>Print (Skriv ut)</b> där diagrammet kan skrivas ut.
Printer Setup (Skrivarinställningar)	Öppnar dialogrutan <b>Printer Setup</b> (Skrivarinställningar) där skrivarinställningarna kan definieras.
Zoom to 100% (Zooma till 100 %)	Ställer tillbaka diagramvisningen till 100 % om zoomfaktorn har ändrats genom att välja rektangulära områden i diagrammet.
Dilution Series Data (Data för spädningsserie)	Visar värdena för skärningspunkter och korrelationskoefficienter.
Properties (Egenskaper)	Välj det här kommandot om du vill ändra egenskaperna för diagramvisningen. Diagrammets titel, axelformat, kurvstil osv., kan anpassas.
Help… (Hjälp)	Öppnar standarddialogrutan hjälp i magellan.

#### 7.4.12 Kontrollfält Kvalitativa resultat

Definiton av cutoff visar en översikt av de aktuellt använda cutoff-intervall och de använda cutoff-gränserna.

I datainformationsfönstret längst ner på skärmen visas en lista över definitioner av cutoffs (försöksgrupper, indata, cutoff-intervall osv.).

Användaren kan se cutoff-resultaten

- Cutoff definition (Definiton av cutoff) Intervall och gränser för intervallen
- Cutoff results (Cutoff-resultat) Cutoff-resultat för varje brunn
- Cutoff statistic (Cutoff-statistik) Statistik över antalet träffar för varje intervall

#### 7.4.13 Kontrollfält Prov-ID:n

Användaren kan visa data från prov-ID-listan.

#### 7.4.14 Kontrollfält Metodlayout

Användaren kan se metodlayouten, dvs. brunnspositioner, layout, replikatinformation, mätparametrar osv.

#### 7.4.15 Kontrollfält QC-validering

Listan över valideringar innehåller försöksgrupperna, indata för varje valideringsgrupp, valideringsformlerna och deras resultat (TRUE eller FALSE).

Om resultatet är TRUE visar det att valideringskriterierna har uppfyllts för denna specifika platta, medan FALSE är ett ogiltigt resultat för hela plattan.

#### 7.4.16 Kontrollfält Miscellaneous (Övrigt)

Felprotokollet loggar alla fel som uppstår under körningen av en metod. Fel kan uppstå under en mätning eller under beräkningen. Gå igenom felprotokollet innan du använder data och resultat.



#### 7.4.17 Box för färgskala

Om värden som ingår i analysplattan presenteras i olika färger, visas dialogrutan **Color scale (Färgskala)**. Färgerna ger en snabb överblick över mätresultaten för de individuella brunnarna. Färgerna, som de används, beror på de inställningar som definieras i färgschemat.

#### 7.4.18 Sammanhangsberoende meny för en brunn

Genom att högerklicka på en brunn i plattlayouten - när du har valt fliken **Edit Method (Redigera metod)** - visas en sammanhangsberoende meny.

#### **Dialogrutan Summary... (Sammanfattning)**

Dialogrutan **Summary (Sammanfattning)** ger en översikt över alla definierade parametrar för en vald brunn.

Information tree (Informationsträd)	Informationsträdet ger en översikt över alla definierade brunnsbaserade parametrar för den valda brunnen. Trädet fylls i enlighet med tillgänglig information, t.ex. om information om prov-ID-listan finns tillgänglig visas prov- ID, pipetteringsstatus osv. Om rådata är tillgängliga kan ytterligare information visas beroende på mätinställningarna och de anslutna instrumenten, t.ex. för kinetiska mätningar visas tidpunkterna.
Knapparna Left (Vänster), Right (Höger), Up (Upp) och Down (Ner)	Riktningsknapparna kan användas för att välja en annan brunn på plattan samtidigt som dialogrutan förblir öppen.
Expand All/ Shrink All (Expandera allt/förminska allt)	Klicka på den här knappen för att visa informationsträdet expanderat till den högsta nivån eller för att minska till den första nivån.
ОК	Stänger dialogrutan Summary (Sammanfattning).

#### **Dialogrutan Details (Detaljer)**

Dialogrutan **Details (Detaljer)** visar resultaten för flera läsningar per brunn för en brunn.

Menyn <b>File (Fil)</b>	Save & Exit (Spara och avsluta) väljs för att spara alla ändringar och stänga dialogrutan.
Menyn <b>Edit (Redigera)</b>	<ul> <li>Menyn Edit (Redigera) innehåller följande kommandon:</li> <li>Klicka på Copy (Kopiera) för att kopiera de enstaka värdena till urklipp, som sedan kan överföras till ett Windowsprogram med hjälp av funktionen klistra in.</li> <li>Välj Recalculate (Beräkna om) för att tvinga fram en ny beräkning av statistiken längst ner i dialogrutan. Det här alternativet är inaktiverat om Automatic Recalculation (Automatisk beräkning) är inställd.</li> </ul>
	<ul> <li>Om Automatic Recalculation (Automatisk beräkning) är inställd, beräknas statistikdata efter varje dölj-/visa-åtgärd.</li> </ul>



Menyn <b>View (Visa)</b>	<ul> <li>Menyn View (Visa) innehåller följande kommandon:</li> <li>Values (Värden) kan väljas för att visa läsningarna som värden.</li> </ul>
	<ul> <li>Anvand det när älternativet för att snabbt och grövt visa flera läsningar per brunn. Graphic (Grafik) kan väljas för att visa de avlästa värdena som färger. Brightness (Ljusstyrka) använder en färg med olika ljusstyrkor för att visa värdena. Pseudo Colors (Pseudofärger) använder olika färger för att visa värdena. En färgskala med ett reglage visas för att ändra intensiteten.</li> </ul>
Menyn <b>Help (Hjälp)</b>	Klicka på <b>Help (Hjälp)</b> för att öppna dialogrutan <b>Help</b> (Hjälp) i <b>magellan</b> .
Klicka på punkterna	När du klickar på <b>reads (läser)</b> kommer det att dölja/visa dem. När en läsning har dolts sätts värdet inom parentes Om färgdisplayen är aktiverad visas läsningen i vitt med en svart cirkulär kant.

#### **Dialogrutan Edit (Redigera)**

Använd det här kommandot för att redigera uppmätta rådata för en vald brunn.

Textfältet Edit raw data (Redigera rådata)	Ett nytt värde för den valda brunnen kan anges. Data visas sedan med symbolen ~ (se kapitel 7.4.6 Särskilda karaktärer).
Knappen <b>Reset</b> (Återställ)	Återställer brunnen till sitt ursprungliga värde.
Kryssrutan <b>Mask (Dölj)</b>	Utesluter värdet av brunnen från beräkningen. Data anges sedan inom parentes. Välj kryssrutan <b>Mask (Dölj)</b> igen för att inkludera värdet igen.
Knapparna Left (Vänster), Right (Höger), Up (Upp) och Down (Ner)	Riktningsknapparna används för att välja en annan brunn på plattan samtidigt som dialogrutan förblir öppen.
ОК	Stänger dialogrutan Edit (Redigera).

#### Dialogrutan Graph: Multilabel (Diagram: Multipla inmärkningar)

Dialogrutan **Graph (Diagram):** Dialogrutan **Multilabel (Multipla inmärkningar)** visar data för de valda brunnarna. Rådata mot brunnsidentiteter visas.

# Sammanhangsberoende meny för diagram med multipla inmärkningar

Genom att högerklicka på diagrammet visas en sammanhangsberoende meny.

Hårkorsmarkör	En hårkorsmarkör visas för att hjälpa till att placera markören på särskilda punkter i diagrammet.
Copy to Clipboard (Kopiera till Urklipp)	Kopierar diagrammet till urklippet som en bitmapp, som kan överföras till alla Windows-program med hjälp av funktionen klistra in.
Save as bitmap (Spara som bitmapp)	Sparar diagrammet som en bitmappfil.



Print… (Skriv ut)	Öppnar dialogrutan <b>Print (Skriv ut)</b> där diagrammet kan skrivas ut.
Printer Setup… (Skrivarinställning ar)	Öppnar dialogrutan <b>Printer Setup</b> (Skrivarinställningar) där skrivarinställningarna kan definieras.
Zoom to 100% (Zooma till 100 %)	Ställer tillbaka diagramvisningen till 100 % om zoomfaktorn har ändrats genom att välja rektangulära områden i diagrammet.
Properties (Egenskaper)	Välj det här kommandot om du vill ändra egenskaperna för diagramvisningen. Diagrammets titel, axelformat, kurvstil osv., kan anpassas.
Help (Hjälp)	Den här knappen öppnar dialogrutan Help (Hjälp) i magellan.

#### Dialogrutan Graph: Kinetics (Diagram: Kinetik)

Dialogrutan **Graph (Diagram): Kinetics (Diagram: Kinetik)** visar diagrammet för kinetiska mätningar av en eller flera valda brunnar. Förklaringen innehåller de beräknade kinetiska parametrarna.

Klicka på punkter för att dölja/visa dem i det kinetiska diagrammet. När en punkt har dolts justeras linjen i diagrammet automatiskt i enlighet med detta och punkten representeras som en transparent symbol i stället för en fylld färgad symbol.

#### Sammanhangsberoende meny för kinetiska diagram

Genom att högerklicka på diagrammet visas en sammanhangsberoende meny.

Hårkorsmarkör	En hårkorsmarkör visas för att hjälpa till att placera markören på särskilda punkter i diagrammet.
Copy to Clipboard (Kopiera till Urklipp)	Kopierar diagrammet till urklippet som en bitmapp, som kan överföras till alla Windows-program med hjälp av funktionen klistra in.
Save as bitmap (Spara som bitmapp)	Sparar diagrammet som en bitmapp.
Print… (Skriv ut)	Öppnar dialogrutan <b>Print (Skriv ut)</b> där diagrammet kan skrivas ut.
Printer Setup (Skrivarinställningar)	Öppnar dialogrutan <b>Printer Setup</b> (Skrivarinställningar) där skrivarinställningarna kan definieras.
Zoom to 100% (Zooma till 100 %)	Ställer tillbaka diagramvisningen till 100 % om zoomfaktorn har ändrats genom att välja rektangulära områden i diagrammet.



#### Visa kinetiska reducerade data

Beräknade kinetiska data (lutningar, start, minima, maxima) kan visualiseras i diagrammet för den kinetiska kurvan eller de kinetiska kurvorna. Lutningar visualiseras som kurvor och starttider, minima och maxima som skärningspunkter. Displayen kan aktiveras/avaktiveras med hjälp av elementet i sammanhangsberoende menyn. **Show Kinetic Reduced Data (Visa kinetiska reducerade data)** i dialogrutan kinetiska diagram. Det aktuella visningsläget sparas och återanvänds för ytterligare visningar.



Temperature curve (Temperaturkurva)	Det här alternativet visar/döljer en kurva som visualiserar temperaturutvecklingen under kinetiska mätningar (om tillgänglig).
Kinetic parameters… (Kinetiska parametrar)	Öppnar dialogrutan <b>Kinetic parameters (Kinetiska</b> <b>parametrar)</b> och ger möjlighet att ändra de kinetiska inställningarna för den eller de valda brunnarna.
Kinetic Data… (Kinetisk data)	Välj det här menyalternativet för att visa en tabell med alla beräknade kinetiska data för de valda brunnarna.
Skalning av Y- axeln	Ett intervall för Y-axeln kan väljas.
Properties (Egenskaper)	Välj den här knappen för att ändra egenskaperna för visning av diagrammet. Diagrammets titel, axelformat, kurvstil osv., kan anpassas.
Hjälp	Öppnar dialogrutan Help (Hjälp) i <b>magellan</b> .





### **EXPERTENS KUNSKAP**

Det är möjligt att ställa in olika kinetiska parametrar för olika brunnar antingen genom att använda den sammanhangsberoende menyn för en brunn i fönstret plattlayout eller genom att använda den sammanhangsberoende menyn i dialogrutan för kinetiska diagram. När du redigerar de kinetiska parametrarna på fliken Edit method (Redigera metod) ställs de kinetiska parametrarna in på samma sätt för alla brunnar.

#### Dialogrutan Graph: Spectra (Diagram: Spektra)

Dialogrutan Graph: Spektra (Diagram: Spektra) visar spektrumet för en skanningsmätning.

#### Sammanhangsberoende meny för spektradiagram i 2D

Genom att högerklicka på diagrammet visas en sammanhangsberoende meny.

Hårkorsmarkör	En hårkorsmarkör visas för att hjälpa till att placera markören på särskilda punkter i diagrammet.
Copy to Clipboard (Kopiera till Urklipp)	Kopierar diagrammet till urklippet som en bitmapp, som kan överföras till alla Windows-program med hjälp av funktionen klistra in.
Save as Bitmap (Spara som bitmapp)	Sparar diagrammet som en bitmapp.
Save as JCAMP-DX (Spara som JCAMP- DX)	Välj det här kommandot för att spara diagrammet som en JCAMP-DX 4.24-fil.
Save as ASCII (Spara som ASCII)	Välj den här knappen för att spara diagrammet som en tabbseparerad ASCII-fil.
Print… (Skriv ut)	Öppnar dialogrutan <b>Print (Skriv ut)</b> där diagrammet kan skrivas ut.
Printer Setup… (Skrivarinställningar)	Öppnar dialogrutan <b>Printer Setup</b> (Skrivarinställningar) där skrivarinställningarna kan definieras.
Zoom to 100% (Zooma till 100 %)	Ställer tillbaka diagramvisningen till 100 % om zoomfaktorn har ändrats genom att välja rektangulära områden i diagrammet.
Spectra Data Reduction (Reducering av spektradata)	Välj det här kommandot för att ändra parameterinställningar för <b>Spectra Data Reduction</b> (Reducering av spektradata).
Spectra Reduced Data (Spektrareducerad data)	Välj det här kommandot för att visa en lista över reducerade spektradata. <b>Anmärkning!</b> Detta kommando aktiveras endast när det finns reducerade spektradata av typvärde tillgängligt.
Data as Table (Data som tabell)	Öppnar en dialogruta som visar spektrumet som en tabell. Detta gör det möjligt att kopiera data till Excel.

Spectrum masked (Spektrum dolt)	Spektrumet och de reducerade data visas som dolda. Detta förhindrar ytterligare beräkningar av denna brunn. Alternativet är endast tillgängligt när råspektrum visas.
Show Prim. Spectrum (Visa prim. spektrum)	Det valda spektrumet och råspektrumet visas tillsammans. Alternativet är endast tillgängligt när reducerat spektrum visas.
Properties (Egenskaper)	Välj det här kommandot om du vill ändra egenskaperna för diagramvisningen. Diagrammets titel, axelformat, kurvstil osv., kan anpassas.
Help (Hjälp)	Öppnar dialogrutan Help (Hjälp) i magellan.



Anmärkning
För att exportera skanningsresultat går du till den
sammanhangsberoende menyn genom att högerklicka på
spektradiagrammet. Välj Data as Table (Data som tabell) i dialogrutan
och kopiera data till Excel eller spara dem som ASCII-fil.

# 7.5 Fliken Edit Method (Redigera metod)

Klicka på den här fliken för att byta till den aktuella metoden och dess inställningar. Varje ändring av metoden resulterar i en ny beräkning av alla data när du går tillbaka till fliken Utvärdera resultat. Dessa ändringar kan sparas i arbetsytan, men de kommer inte att tillämpas på den ursprungliga metodfilen. Se kapitel 4 Create/Edit a Method Wizard (Guiden Skapa/redigera en metod) för mer information

Fliken Edit method (Redigera metod) visas endast om den aktuella användaren har lämpliga rättigheter (kapitel 9.6 Användarrättigheter).



# 7.6 Spara de utvärderade resultaten

Klicka på Next (Nästa) i dialogrutan Results (Resultat) för att nå dialogrutan Save in (Spara i)

Grupprutan <b>Save in</b> (Spara i)	Textfältet <b>Filename</b> (Filnamn): ett standardnamn för filen visas, men kan ändras om så önskas.
	Textfältet <b>File remarks</b> (Filkommentarer): ange kommentarer vid behov.
Grupprutan Audit trail comment (Audit trail kommentar)	Textfältet <b>Audit trail comment</b> (Audit trail kommentar): ange kommentarer till audit trail vid behov (endast tillgängligt för <b>magellan</b> Tracker).
Knappen <b>Save</b> (Spara)	Sparar arbetsytan
<b>Export</b> (Exportera)	Klicka på den här knappen för att öppna dialogrutan <b>Export</b> (Exportera)
Knappen <b>Audit</b> Trail:	Används för att visa <b>Audit trail</b> av filen (endast tillgänglig för <b>magellan</b> Tracker). Se kapitel 4.4 Spara metoden - Audit trail för mer information.
Knappen <b>Signatures</b> (Underskrifter)	Används för att visa de underskrifter som är bifogade till denna fil (endast tillgängligt för <b>magellan</b> Tracker). Se kapitel 4.4 Spara metoden för mer information.
Sign this workspace now (Signera denna arbetsytan nu)	Välj <b>Sign this workspace now</b> (Signera denna arbetsytan nu) för att öppna guiden <b>Attach Signature</b> (Bifoga underskrift) (endast tillgänglig för <b>magellan</b> Tracker).

Klicka på **FINISH (Slutför)** för att stänga guiden **Evaluate Results (Utvärdera resultat).** Om ändringarna inte har sparats ombeds användaren att spara dem.

Klicka på Save (Spara) för att spara data utan att stänga guiden.
## 8. Guiden Bifoga underskrift

## 8.1 Inledning

Flera underskrifter kan appliceras på en post och de finns alltid med i den utskrivna rapporten. Signerade poster kan endast ändras av användare med lämpliga rättigheter. Det är möjligt att helt kontrollera användningen av metoder genom att tillåta användare att endast köra signerade metoder. Endast **Method** (Metod) och Workspace (Arbetsyta) kan signeras.



Anmärkning Den här guiden är endast tillgänglig med magellan Tracker.

Klicka på Attach signature (Bifoga underskrift) för att starta guiden Attach signature (Bifoga underskrift).

Efter dialogrutan Welcome (Välkommen), visas dialogrutan **Select a File (Välj en fil)**. Välj en **Method (Metod)** eller **Workspace (Arbetsyta)** för att signera.

Kombinationsrutan **Show** (Visa)

I kombinationsrutan **Show** (Visa) kan den visade listan över filer ändras i enlighet med urvalet. Möjliga val är:

- All files (Alla filer)
- Unsigned files (Osignerade filer)
- Signed files (Signerade filer)

## 8.2 Signera en fil

Klicka på Next (Nästa) och fönstret Sign (Signera) visas.

Sign		
Signature		
File:	ELISA	
Meaning:	Review     Approval     Custom:	
Comment:	for test purposes	
User Name:	Admin	
Password:	•••••	
Help Cancel	< Back	SIGN
		© 2004 Tecan



<b>Review</b> (Granskning)	Kan endast signeras av en användare med lämpliga rättigheter.
Approval (Godkännande)	Kan endast tillämpas av en användare med lämpliga rättigheter för poster med en redan befintlig granskningsunderskrift. Granskning och godkännande kan inte signeras av samma användare. Inga ändringar får göras i filen mellan granskning och godkännande.
Custom (anpassad)	Ange en anpassad underskriftsbetydelse i textfältet.

Standardanvändarinställningarna tillåter endast administratörer att bifoga underskrifter till granskningar och godkännanden och att ändra signerade filer.

Kommentarer kan läggas till i textfältet Comment (Kommentarer).

I textfältet **User Name (Användarnamn)** måste användarnamnet för den inloggade användaren anges. I textfältet **Password (Lösenord)** måste lösenordet för den inloggade användaren anges.

Klicka på **Finish (Slutför)** för att bekräfta den angivna informationen och signera posten.



#### Anmärkning

Beroende på standardrutinerna för det företag som använder denna programvara kan denna underskrift betraktas som juridiskt bindande. Därför är det mycket viktigt att användarna håller sina lösenord hemliga.

# 9. Ikonen Miscellaneous (Övrigt)

Klicka på ikonen **Miscellaneous (Övrigt)** i guiden och välj mellan följande åtgärder/definitioner:



## 9.1 Instrument control (Instrumentkontroll)

Se avsnitt 3Instrumentkontroll och inställningar..



## 9.2 File handling (Filhantering)

## 9.2.1 Arkivfiler

Grupprutan **Archive Files (Arkivfiler)** ger en översikt och fullständig kontroll över säkerhetskopiering av filer. Den listar alla filer (arbetsytor, metoder, prov-ID-listor eller tillfälliga filer) som sparats av **magellan**.

📶 File handling				
Archive files				
Archive path:				
		(		Archive
File Name	Date:	Remarks	~	Select All
10012008-003.wsp	1/17/2008 12:50:06 PM 1/17/2008 12:55:26 PM	1,25 Diydroxy Vitamin D		Refresh
Vitamin D_3.wsp	1/17/2008 12:55:26 PM			Filter
Vitamin D_1.wsp	1/17/2008 1:20:29 PM			Current Filter:
10012008-003_2.wsp	1/17/2008 1:51:33 PM 1/17/2008 1:51:33 PM			*,*
Elisa_1.mth Elisa_2.mth	1/18/2008 1:03:57 PM 1/18/2008 1:03:57 PM			
() 18012008-001.wsp () 18012008-002.wsp	1/18/2008 1:04:33 PM 1/18/2008 1:04:33 PM		=	
InterfaceTest-001#1	1/18/2008 2:27:11 PM	Endpoint measurement Automation check		
15012008-004_Dain	1/21/2008 1:32:17 PM	Hatomadorrendex		
15012008-004_Dain	1/21/2008 2:06:46 PM		~	
0 of 28 selected		<u>(</u>		
Convert				
Convert from	Convert to			
Import	Logfile			
Import raw data	Save Logfiles			Close

De filer som ska säkerhetskopieras måste väljas från listan **File Name (Filnamn).** Klicka på **Archive (Arkiv)** för att flytta alla valda filer till den angivna säkerhetskopieringskatalogen i lämpliga undermappar.

När filer har attributet **read-only (skrivskyddad)** visas en varningsruta innan filerna raderas.

Grupprutan Archive Files (Arkivfiler) innehåller följande element:

Archive path (Arkivsökväg)	Definierar arkivsökvägen. Klicka på knappen Bläddra 🛄 för att ändra sökvägen.
Listan <b>File Name</b> (Filnamn)	• Listar alla filer som sparats av <b>magellan</b> , filtrerade enligt de alternativ som anges i knappen <b>Filter</b> .
Knappen <b>Archive</b> (Arkiv)	Klicka på <b>Archive</b> (Arkiv) för att flytta alla filer som är markerade i listan <b>File name</b> (Filnamn) till säkerhetskopian (funktion för att klippa och klistra in).

Knappen <b>Select All</b> (Välj alla)	Klicka på <b>Select All</b> (Välj alla) för att välja alla filer som visas i listan <b>File name</b> (Filnamn).
Knappen <b>Refresh</b> (Uppdatera)	Klicka på <b>Refresh</b> (Uppdatera) för att avbryta den aktuella <b>Filter</b> -alternativet och uppdatera fillistan.
Knappen <b>Filter</b>	Den här knappen används för att endast visa specifika filer eller filer med en viss sträng i filnamnet:
	• Ett frågetecken ? kan användas som wildcard för ett enskilt tecken.
	<ul> <li>En asterisk * kan användas som wildcard för inga eller många tecken.</li> </ul>

#### Automatic Archiving (Automatisk arkivering)

Funktionen automatisk arkivering kopierar arbetsytor, metod- och provfiler, som antingen sparas automatiskt eller manuellt av användaren, till en användardefinierad katalog.

Alternativet Automatic Archiving (Automatisk arkivering) är endast tillgängligt i magellan Tracker och kan endast aktiveras av användare med rättigheterna Modify general options (Ändra allmänna alternativ).

ptions				×
Paths Cop	y Plate View	Miscellaneous		
Wizard	welcome page	Start at star	tup I list dialog box tes dialog box	
Demo Move Supp	mode allowed plate in after sta ress 'not monoto	<b>rtup</b> ne' warning for stan	dard curve	
Set next o	counter value Inter: 1	Counter with daily	yreset: 11	
Language	•			
Select:	English		Set	
Additiona	d environments ally accept files fr	om PCs (HUID):	Add HUIDs	
Automatic	archiving le archiving			
Change	password	Options and defau	lt settings audit trail	
		OK Car	ncel Help	
uppruta utomatio	n c	<ul> <li>Kryssruta Markera arkivering</li> </ul>	an <b>Enable arc</b> denna kryssru g.	chiving (Aktivera arkiver uta för att aktivera autom

archivingarkivering.(Automatisk<br/>arkivering)• Klicka på knappen för filbläddraren ("...") för att söka<br/>efter arkivkatalogen.



Anmärkning Det kan hända att en vald arkivkatalog blir ogiltig, t.ex. en

nätverksenhet. I ett sådant fall varnar magellan inte uttryckligen användaren, men gör en notering i loggfilen för systemets Audit Trail.

TECAN.



Anmärkning

Observera att filerna sparas i arkivkatalogen först från och med det ögonblick då funktionen Automatisk arkivering aktiveras. Filer som har sparats innan funktionen Auto Archive (Automatisk arkivering) aktiverades sparas inte i arkiveringskatalogen.

#### 9.2.2 Importera rådata

Genom att använda funktionen **ASCII File Import (Import av ASCII-filer)** kan användaren behandla ASCII-datafiler i **magellan** som redan innehåller mätdata och ytterligare transformationer. Vissa dataformatdefinitioner måste göras i förväg för att garantera korrekt datahämtning. Till exempel måste individuella data skiljas från datalistan med angivna avskiljare.

Ytterligare alternativ som måste specificeras är följande: Plattformat, antal mätningar, mätintervall och typ av dataordning, orientering och fysisk enhet för värdena.

Dataimporten misslyckas om ASCII-filen inte innehåller så många poster som anges i plattformatet.

När du har definierat de nödvändiga alternativen väljer du den önskade filen från fältet **files selection (Urval av filer)** och hämta filinformationen genom att klicka på knappen **Open (Öppna)**.

Dialogrutan **Open (Öppna)** är en standard Windows-dialogruta för att öppna filer med filnavigeringselement, ett textfält för att ange ett filnamn och en rullgardinsmeny för att välja en filtyp (\*.asc).

Dessutom innehåller dialogrutan **Öppna** följande element för att ställa in ASCIIalternativ:

Alternativknappen <b>Format</b>	Den definierar det förväntade dataformatet <b>Table</b> (Tabell) eller <b>Matrix</b> (Matris)
Grupprutan <b>Separator</b> (Avskiljare)	<ul> <li>Den definierar vilket tecken som används för att separera datafälten i listan eller matrisen:</li> <li>Tabulator om <b>Tabulator</b> är valt.</li> <li>Alla andra tecken som anges i textfältet <b>Other</b> (Annat).</li> </ul>
Alternativknappen <b>Orientation</b> (Orientering)	Den definierar orienteringen av data, dvs. antingen horisontell eller vertikal.
Measurement Parame	ters (Mätparameterar)
Info	l rutan visas följande parametrar:

1	rutan visas följande parametrar:
•	Plate format (Plattformat)
•	Data unit (Dataenhet)
•	Number of measurements (Antal mätningar)
•	Measurement interval (Mätningsintervall)



Knappen Measurement **Parameters** (Mätparameterar) Välj mättyp, plattyp, antal kinetikcykler och intervall.

#### 9.2.3 Konvertera till

Dialogrutan Convert Documents - Convert to (Konvertera dokument -Konvertera till) möjliggör användaren att konvertera magellan-filer från den aktuella magellan-versionen till filer för tidigare magellan-versioner.



Anmärkning Det här alternativet är endast tillgängligt om den aktuella användaren har lokala administratörsrättigheter för Windows-systemet.



## FÖRSIKTIGHET

EFTER KONVERTERINGEN MÅSTE METODERNA VALIDERAS, EFTERSOM INNEHÅLLET I DE KONVERTERADE FILERNA KAN SKILJA SIG NÅGOT FRÅN DEN URSPRUNGLIGA FILEN (T.EX. INNEHÅLLET I DEN UTSKRIVNA RAPPORTEN OSV.)

Dialogrutan Convert Documents (Konvertera dokument) innehåller följande element:

Grupprutan <b>Document type</b> (Dokumenttyp)	Välj den dokumenttyp som ska konverteras
Knappen <b>Select all</b> (Välj alla)	Klicka på den här knappen för att välja alla filer som visas i fillistan
Listan <b>Filename</b> (Filnamn)	Alla filer som är tillgängliga för konvertering visas. Klicka på en fil för att välja den för konvertering. Genom att hålla in Shift-tangenten kan du välja flera filer.
Redigeringsfältet <b>Source path</b> (Källsökväg)	Redigeringsfältet <b>Source path</b> (Källsökväg) innehåller sökvägen där de filer som ska konverteras lagras. Klicka på knappen till höger om textfältet för att ändra sökvägen.
Redigeringsfältet <b>Destination path</b> (Destinationssökvä g)	Redigeringsfältet <b>Destination path</b> (Destinationssökväg) innehåller sökvägen där de konverterade filerna lagras. Klicka på knappen till höger om textfältet för att ändra sökvägen.
Knappen <b>Convert</b> document(s) (Konvertera dokument)	Klicka på den här knappen för att starta konverteringen.
Rullgardinsmenyn Version	Välj från rullgardinsmenyn vilken version av <b>magellan</b> som filerna ska konverteras till.
Progress bar (Förloppsindikator)	Denna stapel visar konverteringens förlopp.





## VARNING

#### FILERNA KAN INNEHÅLLA FUNKTIONER SOM INTE ÄR KOMPATIBLA MED DET VALDA FORMATET. DESSA FUNKTIONER ÄR DÅ INTE TILLGÄNGLIGA.

#### 9.2.4 Konvertera från

Dialogrutan **Convert Documents – Convert from (Konvertera dokument – Konvertera från)** möjliggör konvertering av **magellan**-filer från tidigare versioner till det aktuella filformatet. Standard **magellan**-filer kan också konverteras till **magellan** Tracker-filer.



Anmärkning Filer för konvertering visas endast om magellan är ansluten till det instrument (mät- eller demoläge) för vilket filerna genererades.







Anmärkning

Filer från magellan-versioner tidigare än 4.0 måste konverteras för att kunna användas.

Anmärkning Filer från magellan Standard måste konverteras för att kunna användas med magellan Tracker.

## FÖRSIKTIGHET

EFTER KONVERTERINGEN MÅSTE METODERNA VALIDERAS, EFTERSOM INNEHÅLLET I DE KONVERTERADE FILERNA KAN SKILJA SIG NÅGOT FRÅN DEN URSPRUNGLIGA FILEN (T.EX. INNEHÅLLET I DEN UTSKRIVNA RAPPORTEN OSV.)

Dialogrutan Convert Documents (Konvertera dokument) innehåller två flikar.

Fliken magellan 2.x/magellan 3.x innehåller följande element:

Grupprutan <b>Document type</b> (Dokumenttyp)	Välj den dokumenttyp som ska konverteras, t.ex. en metod.
Customize new method specific options (Anpassa nya metodspecifika alternativ) Separator (Avskiljare)	Flera globala alternativ – i <b>magellan</b> -versioner lägre än V4.0 - är nu metodspecifika. För att ställa in dessa alternativ korrekt i metod- och arbetsytorna använder du en av exportknapparna eller knappen Workspace name (Arbetsytans namn)
Knappen <b>Select all</b> (Välj alla)	Klicka på den här knappen för att välja alla filer som visas i fillistan



Listan <b>File name</b> (Filnamn)	Alla filer som är tillgängliga för konvertering visas. Klicka på en fil för att välja den för konvertering. Genom att hålla in Shift-tangenten kan du välja flera filer. <b>Anmärkning!</b> Endast arbetsytor och metoder som skapats med det aktuellt anslutna instrumentet visas.
Redigeringsfältet <b>Source path</b> (Källsökväg)	Redigeringsfältet <b>Source path</b> (Källsökväg) innehåller sökvägen där de filer som ska konverteras lagras. Klicka på till höger om textfältet för att ändra sökvägen.
Redigeringsfältet <b>Destination path</b> (Destinationssökväg)	Redigeringsfältet <b>Destination path</b> (Destinationssökväg) innehåller sökvägen där de konverterade filerna lagras. Klicka på till höger om textfältet för att ändra sökvägen. <b>magellan</b> Tracker: Denna sökväg kan inte ändras.
Knappen <b>Convert</b> document(s) (Konvertera dokument)	Klicka på den här knappen för att starta konverteringen.
Progress bar (Förloppsindikator)	Denna stapel visar konverteringens förlopp.



## VARNING

#### TESTERNA INNEHÅLLER INGA MÄTPARAMETRAR. NÄR ETT TEST OMVANDLAS TILL EN METOD MÅSTE MÄTPARAMETRARNA DEFINIERAS. KONTROLLERA ATT MÄTPARAMETRARNA PASSAR DET VALDA TESTET.

**magellan** 4.x / 5.x-fliken - endast tillgänglig med **magellan** Tracker - innehåller samma fönsterelement som den tidigare beskrivna fliken, förutom att följande alternativ inte är nödvändiga:

• Document type test (Test av dokumenttyp): Testerna är inte tillgängliga med magellan 4.x / 5.x.

Anpassa nya metod<br/>specifika alternativ: Dessa alternativ är redan anpassade i<br/>  $\mbox{magellan}$  4.x / 5.x-filer.

## 9.2.5 Save Log Files (Spara loggfiler)

Alla loggfiler kan sparas som zip-arkiv genom att klicka på knappen **Save Logfiles... (Spara loggfiler)**. Zip-arkivet kan nu namnges och sparas i en angiven katalog. I händelse av mätnings- eller statusfel under utförandet av **magellan** innehåller detta arkiv brunnsdata, status (t.ex. överflöde, "lamp low") eller beräkningsfel och kan enkelt skickas till din lokala kundtjänst för supportbegäran.



## 9.3 Alternativ

Flera val och justeringar som gäller i alla menyer och guider i **magellan** är definierade i dialogrutan **Options (Alternativ)**. Om **magellans** användaradministration är aktiverad, är flera inställningar beroende av användaren. Dialogrutan **Options (Alternativ)** är uppdelad i individuella flikar:

- Fliken Paths (Sökvägar)
- Fliken Copy (Kopiera)
- Fliken Plate View (Plattvy)
- Fliken Miscellaneous (Övrigt)

De begärda alternativen måste anges på motsvarande flikar. Klicka på **OK** för att spara ändringarna och stänga dialogrutan.

Användarberoende alternativ (varje användare kan ändra alternativen för sin egen bekvämlighet) är:

- Standardsökvägar (endast i magellan Standard)
- Alternativ för kopiering till urklipp och Excel
- Plate View (Plattvy)
- (Skip welcome page) Hoppa över välkomstsidan
- Start with wizard list or favorites page (Börja med guide-listan eller favoritsidan)
- Move plate in after startup (Flytta in plattan efter uppstart)

Följande användarberoende alternativ kan inte ställas in i dialogrutan Options (Alternativ):

- Selected Printer (Vald skrivare)
- Printout orientation. (Orientering av utskriften)
- Printout paper size and source. (Pappersstorlek och källa för utskrift)
- Default identifier types. (Standardidentifieringstyper)
- Jumping direction for sample ID input. (Hoppriktning för inmatning av prov-ID)
- Import raw data settings. (Importera inställningar för rådata)
- Standard för Export options (Exportera alternativ), Export to ASCII file (Exportera till ASCII-fil), Export to Excel (Exportera till Excel) och Workspace Name (Arbetsytans namn).
- Graph dialog box size (Storlek på dialogrutan för diagram)

Följande alternativ kan endast ändras av en användare med rätt behörighet. Dessa alternativ gäller sedan för alla användare:

- Language (Språk)
- Demo mode allowed (Demoläge är tillåtet)
- Suppress 'not monotone' warning for standard curve (Undertrycka varningen "not monoton" för standardkurvan)
- Favorites (Favoriter)
- Accepted HUIDs (Accepterade HUID:er)



#### Anmärkning

l magellan Standard kan datasökvägen ställas in av användaren, i magellan Tracker är datasökvägen allmän inställning som gäller för alla användare.



## 9.3.1 Standarddatasökvägar

Den här fliken definierar standardsökvägar för de olika filtyper som används i **magellan** ska sparas i. Dessa fördefinierade sökvägar ställs in vid installationen av programvaran för **magellan** och kan ändras i motsvarande datafält i fliken **Paths (Sökvägar)** 

Def	ault path	5		
Wo	desnace:			
C:\	Users\P	ublic\Docume	ents\Tecan\Magellan\wsp	
Met	hod:			1 10-20-20-20-20-20-20-20-20-20-20-20-20-20
C:\	Users\P	ublic\Docume	ents\Tecan\Magellan\mth	
San	nple ID lis	st:		
C:\	Users\P	ublic\Docume	ents\Tecan\Magellan\smp	
Εхф	ort:			
C:\	Users\P	ublic\Docume	ents\Tecan\Magellan\asc	
Sta	ndard cu	rve:		
C:\	Users\P	ublic\Docume	ents\Tecan\Magellan\wsp	
Bitr	naps:			
C:\	Users\P	ublic\Docume	ents\Tecan\Magellan	
			Change default da	ta path

Fliken Paths (Sökvägar) innehåller följande element:

Textfälten Default paths (Standardsökvägar)	<ul> <li>Ett textfält för var och en av följande filtyper visas med motsvarande standardsökvägar:</li> <li>Workspace (Arbetsyta)</li> <li>Method (Metod)</li> <li>Sample ID list (Prov-ID-lista)</li> </ul>
	• Export (Exportera): för mätdata som exporteras till ASCII-filer.
	<ul> <li>Standard curve (Standardkurva): Sökvägen för standardkurvor är alltid densamma som för arbetsytor.</li> </ul>
	• <b>Bitmaps</b> (Bitmappar): för bitmappar som skapas från dialogrutor för diagram.
🛄 knapparna	Klicka på knapparna bredvid textfälten för att öppna en standard Windows-dialogruta där en annan sökväg för de skapade filerna kan väljas.



Knappen Change default data path (Ändra standarddatasökväg) Finns endast i **magellan** Tracker. Klicka på den här knappen för att ändra standarddatasökvägen för att spara arbetsytor, metoder, prov-ID-listor och standardkurvor. Det krävs rättigheter som administratör och systemadministratör i **magellan** för att ändra standarddatasökvägar.

#### VARNING

PROGRAMVARAN KANSKE INTE KAN TILLDELA RÄTT FILSYSTEMRÄTTIGHETER FÖR EN ANVÄNDARDEFINIERAD STANDARDDATASÖKVÄG! DETTA MÅSTE GÖRAS MANUELLT AV SYSTEMADMINISTRATÖREN. DET FÅR INTE VARA TILLÅTET ATT RADERA ELLER ÄNDRA FILER I STANDARDKATALOGERNA.



Anmärkning I magellan Tracker är det bara export- och bitmappsökningarna som kan ändras.

#### 9.3.2 Alternativ för kopiering/export

I en arbetsyta kan kommandon Edit – Copy (Redigera – Kopiera) och Edit – Copy to Excel (Redigera – Kopiera till Excel) tillåta användaren att exportera mätdata till urklippet eller direkt till ett Excel-kalkylblad. I fliken Copy (Kopiera) anges på vilket sätt innehållet i en platta ska hanteras vid dataöverföring.

Mat	rix format	i		
0	Rows in I	horizontal orde	er / Spectra as rows	
0	Rows in '	vertical order /	/ Spectra as columns	
	Add row/	column labels	(XFluor style)	
Dec	imal cha	racter		
	Point (x.x	x)		
0	Comma (	ktok)		
Con	v to Exc	el ontions		
()	New wor	kbook		
0	New wor	ksheet		
O	nsert into	o worksheet a	t cell A1	
0	Append t	o current work	ksheet	





Fliken Copy (Kopiera) innehåller följande element:

Grupprutan Matrix format (Matrisformat)	<ul> <li>Användaren har också följande möjligheter att definiera hur användaren vill att data ska visas.</li> <li>Alternativknappen Rader i horisontell ordning Raderna med data från mikroplattan kopieras i horisontell ordning.</li> <li>Alternativknappen Rader i vertikal ordning Raderna med data från mikroplattan kopieras i vertikal ordning. Detta resulterar i en överförd matris.</li> <li>Om kryssrutan Add row/column labels (Lägg till etiketter för rad/kolumn) är vald, läggs rad- och kolumninformation som i XFluor till när data kopieras.</li> </ul>
Copy to clipboard options (Alternativ för att kopiera till urklipp)	<ul> <li>överförda data:</li> <li>Alternativknappen Point (x.xxx) (Punkt) Exempel: 7.893</li> <li>Alternativknappen Comma (x,xxx) (Komma) Exempel: 7.893</li> </ul>
Grupprutan Copy to Excel options (Alternativ för att kopiera till Excel)	<ul> <li>Alternativknapparna definierar var överförda data ska placeras i Excel.</li> <li>Alternativknappen New workbook (Ny arbetsbok) Överförda data lagras i det första kalkylbladet i en ny arbetsbok. En ny fil skapas alltid, oavsett om Excel startas eller inte.</li> <li>Alternativknappen New worksheet (Nytt kalkylblad) De överförda data placeras i ett nytt kalkylblad i en öppen, aktiv Excel-arbetsbok. Om Excel inte är öppet när du väljer det här alternativet öppnas det och en ny arbetsbok skapas, och de begärda data infogas i det första kalkylbladet.</li> <li>Alternativknappen Insert into worksheet at cell (Infoga i kalkylbladet i cell) och textfält för cellkoordinater Överförda data placeras i den angivna cellen (standard är cell A1) i ett öppet, aktivt Excel-kalkylblad.</li> <li>Om Excel inte är öppet när du väljer det här alternativet öppnas det och en ny arbetsbok skapas, och de begärda data infogas i det första kalkylbladet.</li> <li>Alternativknappen Insert into worksheet at cell (Infoga i kalkylbladet i cell) och textfält för cellkoordinater Överförda data placeras i den angivna cellen (standard är cell A1) i ett öppet, aktivt Excelkalkylblad.</li> <li>Om Excel inte är öppet när du väljer det här alternativet öppnas det och en ny arbetsbok skapas, och de begärda data infogas i det första kalkylbladet.</li> <li>Alternativknappen Append to current worksheet (Bifoga till aktuellt kalkylblad) Det här alternativet lägger helt enkelt till data i det aktuella kalkylbladet.</li> <li>Om Excel inte är öppet när du väljer det här alternativet öppnas det och en ny arbetsbok skapas, och de begärda data infogas i det första kalkylbladet.</li> </ul>



## 9.3.3 Inställningar för plattvy

I fliken **Plate View (Plattvy)** kan plattvyns utseende anpassas. Färgerna kan ändras för de olika regionerna genom att välja önskad färg från en rullgardinsmeny. Här kan även urvalsläget för märkning av brunnar definieras. Du kan också justera teckensnittsstorleken.

	0	1	2	3	4	
	A					
	В					
	С					Γ
<	-					<b>~</b> ~
Backgrou	nd		Light s	hadow		- -
Plate bott	om	\*  +	Dark s Selecti	hadow ion mode		
Font heigh	nt: med	ium ~			<ul> <li>inters</li> </ul>	ect

Fliken Plate View (Plattvy) innehåller följande element:

Knappen <b>Set</b> (Ställ in)	Bakgrundsfärgen för fönstret <i>Plate view (Plattvy)</i> kan väljas.
Knappen Plate bottom (Plattbotten)	Du kan välja ramfärg för den valda plattan.
Knappen Plate surface (Plattyta)	Ytfärgen på den visade plattan kan väljas.
Rullgardinsmenyn Font height (Teckensnittets höjd)	<ul> <li>Du kan välja teckensnittshöjden för texten i brunnarna:</li> <li>Small (Liten)</li> <li>Medium (Medel):</li> <li>Large (Stor)</li> </ul>
Knappen Light shadow (Ljusskugga)	Du kan välja färgen för ljusskuggan på plattans ram.
Knappen Dark shadow (Mörk kugga)	Du kan välja färgen för mörka skuggan på plattans ram.



Alternativknapparna Selection mode (Urvalsläge) Här fastställs urvalsläget för att markera brunnar med musen: • Union gör det möjligt att välja brunnar endast när

- **Union** gör det möjligt att välja brunnar endast när musen har markerat hela brunnens område.
- Intersect (Skär) gör det möjligt att välja brunnar när endast en del av brunnen har markerats med musen.

## 9.3.4 Övrigt

ptions					×
Paths	Сору	Plate View	Miscellaneous		
Wiz	ard Skip we	lcome page	Start at s Wiz Fav	startup zard list dialog vorites dialog b	box ox
Set Tot	Demom Movepl Suppres next cou	ode allowed ate in after sta s 'not monotor inter value er: 1	itup ne' warning for s Counter with d	tandard curve dailv reset: 1	
Lan Sel	guage lect: E	nglish		~	Set
Reg Ado	gulated e ditionally	nvironments accept files fr	om PCs (HUID):	Add H	UIDs
Auto	omatic a Enable a	rchiving archiving			
Ch	ange pa	ssword	Options and de	fault settings a	udit trail
			ок	Cancel	Help

Fliken Miscellaneous (Övrigt) innehåller följande element:

Grupprutan <b>Wizard</b> (Guide)	När du väljer kryssrutan <b>Skip welcome page (Hoppa</b> över välkomstsidan) visas välkomstsidan för guiden inte när du öppnar en guide. Guiden startar i det andra fönstret.
Grupprutan Start at startup (Starta vid uppstart)	Att välja dialogrutan <b>Favorites (Favoriter)</b> definierar att i stället för <i>Wizard list (Guide-lista)</i> kommer <i>Favorite dialog</i> ( <i>Favoritdialogrutan</i> ) att visas efter start av <b>magellan</b> - programvaran.



Grupprutan	<ul> <li>Kryssrutan Demo mode allowed (Demoläge tillåtet): Välj det här alternativet för att tillåta anslutning till ett simulerat instrument. Detta alternativ är användbart om redan uppmätta data ska utvärderas. Då behövs inget instrument.</li> </ul>
	<ul> <li>Kryssrutan Move plate in after startup (Flytta in plattan efter uppstart): Om kryssrutan är markerad och ett instrument är anslutet flyttas platthållaren in automatiskt efter uppstart.</li> </ul>
	<ul> <li>Kryssrutan Suppress 'not monotone' warning for standard curve (Undertrycka varningen "not monotone" för standardkurvan) Välj det här alternativet om du vill att varningen "not monotone" inte ska visas. Detta alternativ är användbart om flera punkter i en standardkurva tas i ett platt område och det är mycket troligt att monotonin inte är given (dvs. den platta delen av en logaritmisk kurva eller annan kurva där en gräns närmar sig).</li> <li>Anmärkning! Detta alternativ är bara tillgängligt i magellan Tracker.</li> </ul>
Grupprutan	Två alternativ föreslås:
Set next counter value (Ställ in nästa räknarvärde)	<ul> <li>Total counter (Totalräknare) med hjälp av spinnkontrollen kan du ställa in räknaren på valfritt värde.</li> <li>Visar det nummer som kommer att visas i namnet på nästa arbetsyta som skapas.</li> </ul>
	<ul> <li>Counter with daily reset (Räknare med daglig återställning) med hjälp av spinnkontrollen kan du ställa in räknaren på valfritt värde. Visar det aktuella antalet arbetsytor som har skapats under en dag.</li> </ul>
Grupprutan <b>Language (Språk)</b>	Textfältet visar det språk som används. Ett annat språk kan väljas från rullgardinsmenyn. Se <b>Set (Ställ in)</b> nedan.
Knappen <b>Set (Ställ in)</b>	Välj ett annat språk och klicka på <b>Set (Ställ in)</b> för att starta om programmet med det valda språket
Knappen Change password (Byt lösenord)	Klicka på den här knappen för att öppna en dialogruta där användaren kan ändra sitt lösenord (endast tillgängligt när <b>magellan</b> Användaradministration är aktiverad).
Knappen Add HUIDs (Lägg till HUID:er	En lista med ytterligare HUID:er kan definieras så att användaren kan öppna filer som skapats på en annan dator (endast tillgänglig i <b>magellan</b> Tracker).
Automatic Archiving (Automatisk arkivering)	Välj den här funktionen och ange en användardefinierad katalog för att kopiera alla arbetsytor, metod- och prov-ID- filer som antingen sparas automatiskt eller manuellt av användaren till en användardefinierad arkiveringskatalog.
Knappen Options and default settings	Klicka på den här knappen om du vill visa audit trail för alternativ och standardinställningar. Audit trail kan sparas i en ASCII-fil (endast tillgängligt i <b>magellan</b> Tracker).
audit trail… (Alternativ och standardinställning ar audit trail)	Klicka på <b>Save as file (Spara som fil)</b> för att spara informationen som en .log- eller .txt-fil. Texten kan också kopieras till ett ordbehandlingsprogram för utskrift.



# 9.4 User Administration (Användaradministration) – magellan Tracker

Det finns tre typer av administratörer som nämns i denna publikation:

System Administrator (Systemadministratör) är ansvarig för alla ändringar som görs i datorns operativsystem.

Administrator (Administratören) är ansvarig för användarrättigheter i magellan-programvaran.

magellan Administrator (Administratör) har tillgång till alla funktioner i magellan men kan inte ändra användarrättigheter.

**magellan** erbjuder lösenordsskydd för att förhindra missbruk av programvaran och för att begränsa åtkomsten till delar av programvaran baserat på användarrättigheter.



FÖRSIKTIGHET FÖR ATT FÖRHINDRA MISSBRUK AV ANVÄNDARRÄTTIGHETER OCH FÖRFALSKNING AV DATA REKOMMENDERAS DET ATT ANVÄNDARADMINISTRATÖREN INTE HAR MAGELLAN-RÄTTIGHETER (DVS. INTE TILLHÖR EN MAGELLAN-ANVÄNDARGRUPP).

ANVÄNDARADMINISTRATÖREN BÖR HELST VARA EN DEL AV IT-AVDELNINGEN.

Välj miscellaneous (Övrigt) i guiden och klicka på ikonenUser

administration (Användaradministration) Lialogrutan Manage Users and Roles (Hantera användare och roller) visas:

Administrators Magellan_Administrator Magellan_Application_Specialist Magellan_Operator Users		
Full Name	Roles	
Administrator Tom Example	Administrators, Magellan_Adm Magellan_Application_Special	Add User Modify User for Change Password
	nistrator cation_Specialist ator Full Name Administrator Tom Example	nistrator cation_Specialist ator Full Name Roles Administrator Administrators, Magellan_Adm Tom Example Magellan_Application_Special

Användare och roller kan läggas till och ändras, och användarrättigheter kan definieras. Dialogrutan **Manage Users and Roles (Hantera användare och roller)** innehåller följande element:

**Roles** (Roller) Listan **Roles** (Roller) visar alla befintliga användargrupper (användarnivåer).



Add Role (Lägg till roll)	En ny grupp med motsvarande rättigheter kan skapas.
<b>Modify Role</b> (Ändra roll)	Rättigheter för en befintlig grupp kan ändras
<b>Användare</b> (Användare)	Användarlistan visar alla befintliga användare och vilken grupp de tillhör.
Add User (Lägg till Användare)	En ny användare kan skapas.
<b>Modify User</b> (Ändra användare)	Fullständigt namn, användarnivå (grupp) eller lösenord för en befintlig användare kan ändras.
<b>Lock User</b> (Lås användare)	Ett användarkonto kan aktiveras/avaktiveras. Knappen är endast synlig om ett användarnamn är valt.
Audit Trail	Alla ändringar i databasen för användaradministration (t.ex. skapande av grupper/användare, ändring av rättigheter, ändring av alternativ,) registreras i den audit trail för användaradministration som skrivs av servern för användaradministration.
<b>Options</b> (Alternativ)	Allmänna alternativ för inloggning, lösenord och e-post kan ändras
<b>Summary</b> (Sammanfattning)	En detaljerad beskrivning av alla befintliga användare och användargrupper i klartext

#### 9.4.1 Lägg till/ändra användare (magellan Tracker)

Klicka på Add User (Lägg till användare) i dialogrutan Manage Users and Role (Hantera användare och roller) och dialogrutan Create User (Skapa användare) visas:

🤗 Create User		×
Create User	ar name, password and full name of the user to be	
created.	a name, password and full name of the user to be	
<u>U</u> ser Name:	ToEx	1
Full <u>N</u> ame:	Tom Example	1
Password:	××××××× 2	]
<u>C</u> onfirm Password:	жжжжа	
	<u> </u>	

Dialogrutan Create User (Skapa användare) innehåller följande element:

<b>User Name</b> (Användarnamn)	Ett unikt användarnamn måste anges när en ny användare skapas. Namnet kan inte ändras i efterhand.
<b>Full Name</b> (Fullständigt namn)	Ange användarens fullständiga namn - så som det ska visas Detta kan ändras senare.



PasswordAnge det ursprungliga lösenordet. Lösenordet måste ändras<br/>vid nästa inloggning.

#### Klicka på OK och dialogrutan Modify User (Ändra användare) visas:

	User Name:	ToEx
	Full <u>N</u> ame:	Tom Example
	<u>B</u> oles:	Administrators Magellan_Operator
Delete picture	]	Magellan_Administrator

Dialogrutan Modify User (Ändra användare) innehåller följande element:

<b>Full Name</b> (Fullständigt namn)	Ange användarens fullständiga namn - så som det ska visas			
Roles (Roller)	Välj en av de befintliga användargrupperna. Användaren tilldelas de rättigheter som är kopplade till gruppen.			
<b>Change picture</b> (Ändra bild)	Användarens bild kan läggas till eller ändras.			
<b>Delete picture</b> (Radera bild)	Radera användarens bild.			
Change Password (Ändra lösenord)	Ett lösenord för en ny användare kan definieras. Observera att administratören inte behöver ange det gamla lösenordet.			

Om du vill ändra en användare klickar du på **Modify User (Ändra användare)** i dialogrutan **Manage Users and Roles (Hantera användare och roller)**. I dialogrutan **Modify User (Ändra användare)** fortsätter du enligt beskrivningen ovan.

## 9.4.2 Lägg till/ändra roll

Klicka på Add Role (Lägg till roll) i dialogrutan Manage Users and Roles (Hantera användare och roller) och dialogrutan Create New Role (Skapa ny roll) visas. Definiera ett unikt Role Name (Rollnamn). Observera att detta namn inte kan ändras i efterhand. Stäng dialogrutan Add Role (Lägg till roll) genom att klicka på OK. Öppna dialogrutan Modify Role (Ändra roll):

Update Role Update the role by n	nodifying associations with application rig	hts. The role
name cannot be mo <u>R</u> ole Name:	Users	
Application:	Magellan	~
<u>B</u> ights:	<ul> <li>Create sample id lists</li> <li>Edit sample id lists</li> <li>Sign sample id lists</li> <li>Modify signed sample id lists</li> <li>Create methods</li> <li>Edit methods</li> <li>Sign methods</li> <li>Modify signed methods</li> <li>Run unsigned methods</li> </ul>	

Dialogrutan Modify Role (Ändra roll) innehåller följande element:

<b>Role Name</b> (Rollnamn)	Rollnamnet definieras i dialogrutan Add Role (Lägg till roll).
Application (Program)	Tecan-programvarukomponent, dvs. magellan.
<b>Rights</b> (Rättigheter)	Rollen representerar en specifik användarnivå och motsvarande användarrättigheter kan väljas.



## 9.4.3 Audit trail för användaradministration

-----

Klicka på Audit Trail i dialogrutan Manage Users and Roles (Hantera användare och roller) och dialogrutan Audit trail visas:

C:\Progra C:\Docur C:\Docur 2011 2011 2011	amData\Tecan\` nents and Settin 8-02-10 09-28-28-38 8-02-10 09-28-28-66 8-02-10 09-28-29-70 8-02-10 09-28-29-70	Tecan User Mana gs\All Users\App 80.log 61.log 08.log	agement\v1.1\Au Ilication Data\Te	dit\ can\Tecan User	Manageme
201	8-02-10 09-28-58-2	92.log			
	<u>.</u>				3
Q	S = checksum of I	file is correct	S = checksum of fi	le is incorrect	
This selected file w	= checksum of I vas modified	file is correct	= checksum of fi	le is incorrect	
This selected file w	= checksum of l res modified	file is correct	= checksum of fi	le is incorrect	3
: This selected file w by on	= checksum of I vas modified	Rie is correct 😵	= checksum of fi	le is incarrect	

Alla ändringar i databasen för användaradministration registreras i den audit trail för användarhantering som skrivs av servern för användaradministration.

Om du vill visa en loggfil väljer du den i listan (välj kryssrutan) och klickar på knappen **Print (Skriv ut)**. Filen öppnas via en PDF-läsare och kan sedan skrivas ut. Den gröna bocken markerar filer med korrekt kontrollsumma, det röda korset markerar filer med felaktig kontrollsumma.



Anmärkning Observera att den aktuella loggfilen alltid har ett rött kryss eftersom filen fortfarande ändras så länge programmet är öppet.

#### User Rights (Användarrättigheter)

Det finns tre olika säkerhetsnivåer för användarrättigheter, varav den högsta nivån är **magellan Administrator (Administratör)**, som har tillgång till alla programfunktioner. **Programspecialist** och **Operatör**-nivåerna är alltmer begränsade.

magellan Standard: De standardiserade rättigheterna tilldelas enligt nedan.

**magellan Tracker**: Varje rättighet kan överlåtas eller återkallas av en administratör och som standard överlåts den enligt nedan. Standardrättigheterna är baserade på en typisk laboratoriemiljö.



#### FÖRSIKTIGHET ADMINISTRATÖRER HAR ANSVARET FÖR ATT SE TILL ATT INSTÄLLNINGARNA (OCH EVENTUELLA ÄNDRINGAR) STÄMMER ÖVERENS MED DEN LABORATORIEMILJÖ SOM DEFINIERAS AV STANDARDISERADE TILLVÄGAGÅNGSSÄTT (STANDARD OPERATING PROCEDURES) OCH ATT DE ÖVERENSSTÄMMER MED

GÄLLANDE LAGAR.



Anmärkning Vid klinisk diagnostik måste den operativa myndigheten validera alla metoder för att garantera resultatens säkerhet. En metod anses validerad när den har undertecknats av den operativa myndigheten. Den operativa myndigheten är helt ansvarig för de resultat som uppnås.

## 9.4.4 Alternativ för användaradministration

Klicka på Options (Alternativ) i dialogrutan Manage Users and Roles (Hantera användare och roller) och dialogrutan Edit Options (Redigera alternativ) visas.

#### Alternativ för Login (Inloggning)

dit Optio	ons			×
Login	Password	E-Mail		
Loc	k application	fter the specified minutes o	fidle time:	
	15			
Exit	application a mpts:	d lock account after the sp	ec <mark>ifie</mark> d number of unsuccessful logir	1
	3	-		
			OK C	ancel

Lock application (Lås program)	Låsning av program efter en viss tid (1 min–36500 h). Ett lösenord måste anges för att låsa upp programmet.
Unsuccessful logins (Misslyckade	Efter ett antal misslyckade inloggningar i följd (1–100) låses användarkontot och eventuellt skickas ett varningsmeddelande till systemadministratören.
inloggningar)	Om användarkontot för en användaradministratör har låsts på grund av misslyckade inloggningar förblir kontot låst under en inaktivitetstid på 45 minuter. Efter denna väntetid låses kontot upp och kan användas igen.







#### FÖRSIKTIGHET

OM LÖSENORDET FÖR ANVÄNDARADMINISTRATÖR HAR GLÖMTS BORT KAN INSTÄLLNINGARNA FÖR ANVÄNDARADMINISTRATION INTE ÄNDRAS. I SÅ FALL BÖR DU FÖRST AVINSTALLERA MAGELLAN OCH DÄREFTER GÖRA EN NY INSTALLATION (SE KAPITEL 1.5 INSTALLATIONSPROCEDUR FÖR PROGRAMVARA).



Anmärkning När du avinstallerar magellan kommer inga befintliga magellan-filer att gå förlorade. Alla tidigare filer finns i en säkerhetskopieringsmapp som skapas under den nya installationen.

#### Alternativ för Password (Lösenord)

lit Optic	ons				>
Login	Password	E-Mail			
Mini	imum passwo	rd length:			
	5	-		Allow empty p	assword
Mini	imum require	number of no	n-alphanumeric chara	cters: (except for Admi	nistrator)
	0	<b></b>			
Mini	imum require	d number of nu	meric digits:		
	0	+			
Mini	imum numbei	of password of	hanges before re-use		
	3	-			
Pas	sword expire	s after the spe	ified number of days:		
	90	-			
				OK	Cancel

Lösenord måste innehålla ett visst minsta antal tecken. Lösenord kan definieras så att de innehåller icke-alfanumeriska tecken och/eller numeriska siffror. Lösenordet kan inte återanvändas.

#### Minimum password length (Minsta lösenordslängd):

Ange det antal tecken som krävs.

Minimum required number of non-alphanumeric characters (Minsta antal icke-alfanumeriska tecken som krävs):

Ange det antal icke-alfanumeriska tecken som krävs.

Minimum required number of numeric digits (Minsta antal numeriska siffror som krävs):

Ange önskat antal siffror.

Minimum number of password changes before re-use (Minsta antal ändringar av lösenordet innan det återanvänds):

Ange antalet lösenordsändringar.

Password expires after the specified number if of days (Lösenordet upphör att gälla efter det angivna antalet dagar):



Användarens lösenord upphör att gälla efter ett konfigurerbart antal dagar. Användaren måste då välja ett nytt lösenord. Ange önskat antal dagar (intervall: 1 till 36500 dagar).

#### Allow empty password (Tillåt tomma lösenord):

Markera den här kryssrutan om det är tillåtet att ha ett tomt lösenord för alla användare utom administratören.

#### Alternativ för Email (E-postmeddelande)

ogin Password E-Mail	
☑ Notify via e-mail	
Mail to:	
Send by SMTP     Server:	
Port:	
Security Protocol: O SSL O TLS	None
Return Address:	
Authentication:	
User.	
Password:	

administratör kan underrättas om potentiella säkerhetsattacker (ett användarkonto har låsts på grund av ett antal misslyckade inloggningar).

Ett e-postmeddelande kan skickas till den angivna e-postadressen:

• **SMTP-server**: SMTP-serverns IP-adress ska anges (t.ex. SEUATEXCH01.eu.tecan.net)

Klicka på **Test Mail (Testa e-post)** för att skicka ett test e-postmeddelande för att kontrollera att inställningarna är korrekta.



Anmärkning Meddelandet via e-post kräver en SMTP-server utan autentisering.

#### 9.4.5 Sammanfattning av användaradministration

En detaljerad beskrivning av alla befintliga användare och användargrupper i klartext kan erhållas genom att klicka på **Summary (Sammanfattning)** i dialogrutan **Manage Users and Roles (Hantera användare och roller)**.

Texten kan kopieras till ett ordbehandlingsprogram för utskrift.

# 9.5 User Administration (Användaradministration) - magellan Standard

Programvaran **magellan** skyddar mot missbruk av programvaran eller mot att obehöriga användare får tillgång till programvaran genom att tillhandahålla ett lösenordsskydd.

ι	Jser Administration		
	User name Admin (Admin) APSP (App.Specialist)	Rights Magellan_Administ Magellan_Operator	Add User Modify Disable
	Password protection (on/off)		Login
		OK Cancel	Help

Användarprofiler kan definieras på tre olika säkerhetsnivåer. Den högsta säkerhetsnivån i programvaran är den som gäller för **magellan Administrator** (Administratör). Denna nivå ger tillgång till alla programmets funktioner och möjligheter att redigera data. För Application Specialist (Programspecialist) och Operator (Operatör)-nivå är alternativen alltmer begränsade.

Endast **User Administrator (Användaradministratören)** har möjlighet att aktivera eller inaktivera programvarans lösenordsskydd, definiera nya användare eller radera gamla användare. Om användarna är inloggade under sitt användarnamn ger systemet endast tillgång till alternativ enligt definitionen av användarrättigheterna.

Dialogrutan **User administration (Användaradministration)** innehåller följande element:

Knappen <b>Login</b> (Inloggning)	Dialogrutan <b>Login</b> (Inloggning) visas. En användare kan logga in med sitt användarnamn.
<b>User list</b> (Användarlista)	<ul> <li>Alla användare som aktuellt är konfigurerade med sina User name (Användarnamn) och Rights (Rättigheter) visas i denna lista.</li> </ul>
	Här kan du välja vilken användare som ska redigeras eller raderas.
Knappen <b>Add User</b> (Lägg till användare)	Knappen <b>Add User</b> (Lägg till användare) öppnar dialogrutan <b>User Rights</b> (Användarrättigheter) där du kan ange lösenord, namn och användar-ID för att definiera en ny användare.
Knappen <b>Modify</b> (Ändra)	Knappen <b>Modify</b> (Ändra) öppnar dialogrutan <b>User Rights</b> (Användarrättigheter) för att ändra den valda användaren i användarlistan.
Knappen <b>Disable</b> (Inaktivera)	Knappen <b>Inaktivera</b> inaktiverar den valda användaren i den aktuella användarlistan.



Kryssrutan	Den här kryssrutan gör det möjligt att slå på/av
Password protection	programvarans lösenordsskydd. Om den är aktiverad, kan
Zoom	dialogrutan Login (Inloggning) visas när du startar
	magellan.

#### 9.5.1 Lägg till/ändra användare (magellan Standard)

Den här dialogrutan gör det möjligt att lägga till eller redigera användare och tilldela dem till en användargrupp. För rättigheter för olika användargrupper se kapitel 9.6 Användarrättigheter .

Dialogrutan User Rights (Användarrättigheter) innehåller följande element:

Textfältet <b>User Name</b> (Användarnamn)	Ange användarens fullständiga namn – så som det ska visas
Textfältet <b>UserID</b> (Användar-ID)	Definiera det användar-ID som användaren måste ange för inloggning
Alternativknapparna <b>User Rights</b> (Användarrättigheter)	<ul> <li>Definiera nivån på användarens tillgång till programvarans funktioner:</li> <li>Operator (Operatör)</li> <li>Application Specialist (Programspecialist)</li> <li>Administrator (Administratör)</li> </ul>
Textfältet Password (Lösenord)	Ange användarens lösenord (minst 5 tecken)
Textfältet <b>Confirm password</b> (Bekräfta lösenordet)	Ange lösenordet på nytt. Av säkerhetsskäl visas lösenordstecknen som symboler i det här fältet.

Klicka på knappen **OK** för att spara och stänga dialogrutan.

Klicka på knappen Avbryt för att stänga dialogrutan utan att spara inställningarna.



Om inga användare har definierats alls och användaradministrationen ska användas måste en administratör först definieras. När du klickar på inloggningsknappen visas därför ett kort meddelande. Därefter visas dialogrutan Användarrättigheter för att ange administratörens användarinformation (se kapitel 2.3 Starta magellan/Inloggning).

## 9.5.2 Inloggning

Den här knappen är endast aktiverad om användaradministrationen är aktiverad, men **Password protection (Lösenordsskydd)** inte har valts. Därför visades dialogrutan **Inloggning** inte vid uppstart. För att aktivera den lösenordsskyddade användaradministrationen klickar du på knappen **Login (Inloggning)**.

## 9.5.3 Ändra användare

Om du vill ändra den aktiva användaren klickar du på ikonen Change User

(Ändra användare) i guiden. Dialogrutan Login (Inloggning) visas.





## 9.6 Användarrättigheter

Det finns tre olika säkerhetsnivåer för användarrättigheter, varav den högsta nivån är Administrator (Administratör), som har tillgång till alla programfunktioner. Application Specialist (Programspecialist) och Operator (Operatör)-nivåerna är alltmer begränsade.

magellan Standard: De standardrättigheterna tilldelas enligt nedan.

**magellan Tracker**: Varje rättighet kan tilldelas eller återkallas av en användaradministratör och som standard tilldelas den enligt nedan. Standardrättigheterna är baserade på en typisk laboratoriemiljö.





## FÖRSIKTIGHET

ANVÄNDARADMINISTRATÖRER HAR ANSVARET FÖR ATT SE TILL ATT INSTÄLLNINGARNA (OCH EVENTUELLA ÄNDRINGAR) STÄMMER ÖVERENS MED DEN LABORATORIEMILJÖ SOM DEFINIERAS AV STANDARDISERADE TILLVÄGAGÅNGSSÄTT (STANDARD OPERATING PROCEDURES) OCH ATT DE ÖVERENSSTÄMMER MED GÄLLANDE LAGAR.

#### Anmärkning

Vid klinisk diagnostik måste den operativa myndigheten validera alla metoder för att garantera resultatens säkerhet. En metod anses validerad när den har undertecknats av den operativa myndigheten. Den operativa myndigheten är helt ansvarig för de resultat som uppnås.

Användarrättigheterna sammanfattas i tabellen nedan och gäller för både **magellan** -versioner, dvs. **magellan Tracker** och **magellan Standard**, om inget annat anges:



	magellan Administrator	magellan Application Specialist (Program-	magellan Operator
User Rights (Användarrättigheter)	(Administrator)	specialist)	(Operator)
Sample ID List (Prov-ID-lista)			
Create/Edit a Sample ID List (Skapa prov- ID-listor)	Ja	Ja	Ja
Skapa en ny prov-ID-lista eller spara en befintlig lista med ett nytt namn.			
Edit sample ID lists (Redigera prov-ID-listor)			
Redigera en befintlig prov-ID-lista och spara den med samma namn.	Ja	Ja	Nej
Sign sample ID lists* (Signera prov-ID-listor)			
Signera en befintlig prov-ID-lista för granskning/godkännande.	Reserverad för framtida användning.		rändning.
<b>Modify signed sample ID lists</b> * (Ändra signerade prov-ID-listor)	Reserverad för framtida användning.		röndning
Redigera eller ändra prov-ID:n i en signerad prov-ID-lista.			andning.
Method (Metod)			
Create methods (Skapa metoder)			
Skapa en ny metod eller spara en befintlig metod med ett nytt namn.	Ja	Ja	Nej
Edit methods (Redigera metoder)			
Redigera en befintlig metod och spara den med samma namn.	Ja	Ja	Nej
Sign methods* (Signera metoder)			
Signera en befintlig metod för granskning/godkännande.			
Anmärkning: Rätt <b>Sign methods</b> (Signera metoder) räcker inte för att signera metodfiler. Du behöver också rätt <b>Edit methods</b> (Redigera metoder).	Ja	Nej	Nej
Dessutom behövs rättigheten <b>Modify signed</b> <b>methods</b> (Ändra signerade metoder) för att lägga till <b>Approval signature</b> (Underskrift för godkännande) till filer som redan har en underskrift för <b>Review</b> (Granskning).			
<b>Modify signed methods</b> * (Ändra signerade metoder)	Ja	Nei	Nei
Redigera en signerad metod och spara den med samma namn.		-,	-,



User Rights (Användarrättigheter)	<b>magellan</b> <b>Administrator</b> (Administratör)	magellan Application Specialist (Program- specialist)	<b>magellan</b> <b>Operator</b> (Operatör)
Run unsigned methods* (Kör osignerade metoder) Endast med denna rättighet kan endast godkända metoder användas. Anmärkning: Denna rättighet kan användas i kombination med rättigheten Run not approved methods (Kör inte godkända metoder) för att tillåta användaren att köra alla metoder.	Ja	Ja	Nej
<b>Setup favorites</b> (Ställ in favoriter) Lägg till/ta bort metoder från favoritlistan.	Ja	Nej	Nej
<b>Define multilabel measurement</b> (Definiera mätning med multipla inmärkningar) Definiera metoder med multipla inmärkningar	Ja	Ja	Nej
<b>Define kinetic transformations</b> (Definiera kinetiska transformationer) Definiera metoder som använder kinetiska transformationer.	Ja	Ja	Nej
<b>Define concentration transformations</b> (Definiera koncentrationstransformationer) Definiera metoder som använder koncentrationstransformationer.	Ja	Ja	Nej
<b>Define alias</b> (Definiera alias) Definiera metoder som använder alias för identifierare.	Ja	Ja	Nej
Workspace (Arbetsyta)			
Skapa arbetsytor Utför en mätning och spara arbetsytan med ett nytt namn – eller – spara en befintlig arbetsyta med ett nytt namn.	Ja	Ja	Ja
Edit workspaces (Redigera arbetsytor) Redigera en befintlig arbetsyta och spara det med samma namn.	Ja	Ja	Nej



User Rights (Användarrättigheter)	<b>magellan</b> <b>Administrator</b> (Administratör)	magellan Application Specialist (Program- specialist)	<b>magellan</b> <b>Operator</b> (Operatör)
Sign workspaces* (Signera arbetsytor) Signera en befintlig arbetsyta för granskning/godkännande. Anmärkning: Rättigheten Sign workspaces (Signera arbetsytor) räcker inte för att signera arbetsytor. Du behöver också rättigheterna Edit workspaces (Redigera arbetsytor). Dessutom kan rättigheten Edit signed workspaces (Redigera signerade arbetsytor) behövs också för att lägga till Approval signature (Underskrift för godkännande) till filer som redan har överskriften Review (Granskning).	Ja	Nej	Nej
Edit signed workspaces* (Redigera undertecknade arbetsytor) Redigera en signerad arbetsyta och spara det med samma namn.	Ja	Nej	Nej
<b>Continue evaluation when error</b> (Fortsätt utvärderingen när fel uppstår) Fortsätt beräkningen efter att fel har uppstått	Ja	Ja	Nej <sup>1</sup> /Ja²
<b>Modify raw data (samples)</b> (Ändra rådata - prover) Redigera eller dölj mätdata för prover.	Ja	Nej	Nej
Modify raw data (standards) (Ändra rådata - standarder) Redigera eller dölj mätdata för standarder, kontroller och blanka.	Ja	Nej	Nej
Import raw data (Importera rådata - standarder) Importera rådata från en ASCII-fil i en befintlig arbetsyta.	Ja	Nej	Nej
<b>Modify evaluation</b> (Ändra utvärdering) Ändra inställningar för utvärdering (metod) i en befintlig arbetsyta	Ja	Ja	Nej
Reevaluate with another method (Omvärdera med en annan metod) Omvärdera en befintlig arbetsyta med en annan metod.	Ja	Ja	Nej
Modify layout (Utforma layout) Ändra mätparametrar, koncentration och layout innan du påbörjar en mätning.	Ja	Ja	Nej



User Rights (Användarrättigheter)	<b>magellan</b> <b>Administrator</b> (Administratör)	magellan Application Specialist (Program- specialist)	<b>magellan</b> <b>Operator</b> (Operatör)
<b>Create standard curve files</b> (Skapa filer med standardkurvor) Skapa en ny standardkurva och spara den med ett nytt namn.	Ja	Ja	Nej
Edit standard curve files (Redigera filer med standardkurvor) Redigera en befintlig standardkurva och spara den med samma namn	Ja	Ja	Nej
Sign standard curve files* (Signera filer med standardkurvor) Signera en befintlig standardkurva för granskning/godkännande.	Reserve	rad för framtida anv	rändning.
Modify signed standard curve files* (Ändra signerade filer med standardkurvor*) Redigera och ändra inställningar i en signerad fil med standardkurva.	Reserverad för framtida användning.		
General (Allmänt)			
<b>Archive files</b> (Arkivfiler) Flytta filer till en arkivplats	Ja	Nej	Nej
<b>Set default paths</b> ** (Ange standardsökvägar) Ange individuella sökvägar för att spara arbetsyta, metod, prov-ID-lista, export- och bitmappfiler.	Ja	Ja	Ja
<b>Define filter slides</b> (Definiera filterhållare) Ändra filterdefinitionen för instrumentet	Ja	Nej	Nej
Modify general options (Ändra de allmänna alternativen) Välj språk, undertryck inte monotona varningar, tillåt demoläge.	Ja	Nej	Nej
Print unsigned workspaces* (Skriv ut arbetsytor utan underskrift) Resultaten av en arbetsyta utan underskrift kan skrivas ut.	Ja	Ja	Ja
Run not approved methods (Kör inte godkända metoder) Denna rättighet ger användaren möjlighet att använda redan granskade, men ännu inte godkända metoder och godkända metoder. Anmärkning: Denna rättighet kan användas i kombination med rätten att Run unsigned methods (Köra osignerade metoder) för att tillåta användaren att köra alla metoder.	Ja	Ja	Nej

User Rights (Användarrättigheter)	<b>magellan</b> <b>Administrator</b> (Administratör)	magellan Application Specialist (Program- specialist)	<b>magellan</b> <b>Operator</b> (Operatör)
Edit reviewed method* (Redigera granskad metod) Granskade (men inte godkända) metoder kan redigeras.	Ja	Nej	Nej
Edit reviewed workspace* (Redigera granskad arbetsyta) Granskade (men inte godkända) arbetsytor kan redigeras.	Ja	Nej	Nej
Reader Server (Server för läsare)			
Save plate definition files (Spara plattdefinitionsfiler) Skapa eller redigera plattdefinitionsfiler.	Ja	Nej <sup>1</sup> /Ja <sup>2</sup>	Nej
Save spin profile (Spara spinnprofil) Skapa eller redigera filer för spinnprofiler.	Ja	Nej	Nej

\* användarrättigheter som endast finns tillgängliga i magellan Tracker

\*\* användarrättigheter som endast finns i magellan Standard

<sup>1</sup> magellan Tracker

<sup>2</sup> magellan Standard

## 9.7 Information om magellan

Välj **About magellan (Information om magellan)** för att visa information om upphovsrätt och ansvarsfriskrivning för programvara:

I dialogrutan **About magellan (Information om magellan)** visas information om den aktuellt installerade versionen av **magellan**, komponenternas versionsnummer (fliken **components (Komponenter)**), UDI-information och användarens programlicens (fliken **license (Licens)**).

#### **Register Wizard (Guiden Registrera)**

För ytterligare information om **Register magellan Wizard (Guiden Registrera magellan)** se kapitel 2.4.4 Registreringsguiden.



## 10. Ytterligare funktioner för magellan Tracker

## **10.1** User administration (Användaradministration)

Med **magellan** Tracker är användaradministrationen obligatoriskt. Se kapitel 9.4 User Administration (Användaradministration) – magellan Tracker för mer information.

#### 10.1.1 Audit Trail

För att uppfylla föreskrifterna i **FDA 21 CFR part 11** dokumenteras varje steg som görs i **magellan** Tracker. I **Audit trail** listas alla ändringar som rör följande data:

- en metodfil
- en arbetsyta
- en prov-ID-lista
- en fil med en standardkurva
- databasen för användaradministration eller
- alternativ och standardinställningar.

Varje inmatning består av användaren (namn och fullständigt namn), datum och tid för ändringen, om filen skapades eller ändrades och eventuell kommentarer till audit trail.

#### Audit trail för metoder, arbetsytor och prov-ID-listor

Du kan visa audit trail med hjälp av knappen **Audit trail** på sidan Save (Spara) i guiderna.

#### Audit trail av en standardkurva

Standardkurvor kan sparas och sedan användas för utvärdering av data från en annan arbetsyta. Funktionen för att ladda en standardkurva från en extern fil finns i kapitel 4.3.8 Koncentrationer:Standardkurva i guiden Skapa/redigera en metod. För att öppna en tidigare sparad standardkurva i diagrammet för en standardkurva se kapitel 7.4.11 Kontrollfält Koncentrationer:

Om du vill visa audit trail för den laddade standardkurvan väljer du **Audit trail** från menyn **View (Visa)** i dialogrutan för diagrammet för en standardkurva.

#### Audit trail av databasen för användaradministration

Om du vill visa Audit trail för databasen för användaradministration väljer du ikonen **Miscellaneous (Övrigt)** i **Wizard List (Guide-lista)** och klickar på **User Administration (Användaradministration)**.

#### Audit trail av alternativ och standardinställningar

Om du vill visa Audit trail för alla alternativ och standardinställningar väljer du ikonen **Miscellaneous (Övrigt)** i startsidan för **Wizard List (Guide-lista)** och klickar på **Options (Alternativ)**. I fliken **Miscellaneous (Övrigt)** kan knappen **Audit trail** sedan väljas.

## 10.2 Filhantering

De uppmätta data sparas alltid direkt efter mätningen.



Observera att detta endast gäller för magellan Tracker: För att garantera att inga data kan manipuleras (raderas) bör magellananvändare inte vara Windows-administratörer.

TECAN.

## 10.2.1 Spara en fil

En fil som skapas och sparas i **magellan** Tracker, anges med ett "#"-tecken i filnamnet. Det tillagda "#"-tecknet kan endast ses i den lämpliga underkatalogen i **magellan**.

Exempel:

method#1.mth workspace#1.wsp sampleIDlist#1.smp

## 10.2.2 Ändra en metodfil eller en arbetsyta

Varje gång en metodfil eller arbetsyta ändras och sparas ökar numret bakom "#" i filnamnet automatiskt (t.ex.: första versionen av en arbetsyta: name#1.wsp; redigera och spara den nya versionen för att få name#2.wsp; efter nästa ändring skapas name#3.wsp...). Alla versioner av en fil listas i lämpliga underkataloger i **magellan**.



Anmärkning När du överför Tracker-filer måste du överföra alla versioner av filen.

## 10.2.3 Öppna en fil

Endast den senast sparade versionen av en fil är tillgänglig i guiderna för att skapa/redigera en metod, skapa/redigera en prov-ID-lista, utvärdera resultat och starta mätning.

Om du vill visa en tidigare version av en fil öppnar du **Audit trail** på sidan där filen sparas, välj den önskade versionen av filen genom att klicka en gång på numret och klicka på knappen **Save previous version as... (Spara tidigare version som)**. Byt namn på filen och lägg till några filkommentarer eller kommentarer till Audit trail och öppna filen med guiden **Create/Edit a Method (Skapa/redigera en metod)**.



Anmärkning Om du vill jämföra en metod med tidigare versioner måste du göra en utskrift, eftersom två fönster för förhandsgranskning av utskrift inte kan öppnas samtidigt.

## 10.2.4 Öppna en fil som skapats på en annan PC – Lägg till HUID:er

I licensproceduren för magellan finns tre viktiga nummer:

- Programvarans serienummer
- HUID (Hardware Unit Identification Number (Identifikationsnummer för maskinvara)) och
- Licensnummer (se kapitel 2.4.3 Licensiera magellan).



HUID för en persondator konfigureras under registrering av **magellan** och alla filer som skapas och sparas på denna PC kopplas till detta speciella HUID.

Om det finns behov av att arbeta med **magellan** Tracker-filer som skapats på en annan PC måste HUID för den andra PC:n accepteras.

När en användare med rättigheten att **Modify general options (Ändra allmänna alternativ)** försöker öppna ett dokument (prov-ID-lista, metod, arbetsyta) från en annan PC, dvs. med ett annat HUID-nummer, ombeds användaren automatiskt att lägga till det numret i listan över kända HUID-nummer. Om knappen **Accept HUID (Acceptera HUID)** klickas, läggs HUID-numret till i listan och dokumentet öppnas utan ytterligare interaktioner.

#### Detta HUID-nummer accepteras därför alltid.

För att visa listan över accepterade HUID-nummer för andra PC:n klickar du på ikonen **Miscellaneous (Övrigt)** på guide-listans startsida → **Options (Alternativ)** → Fliken **Miscellaneous (Övrigt)**. Klicka på knappen **Add HUID (Lägg till HUID)** och dialogrutan med de andra datorernas HUID-nummer visas.

Invalid HUID	×
This file was created on another computer (HUID: 9408G-3U335FB)! (C:\Users\Public\Documents\Tecan\Magellan\mth\Application Example_en.mth	)
Press 'Accept HUID' to add the given HUID to the list of available HUIDs. Files fr without getting any warning message.	om computers in this list can be opend
	Accept HUID Reject HUID

## 10.3 Systemets Audit Trail

Systemets Audit Trail är endast tillgängligt i **magellan Tracker**-versionen. Systemets Audit Trail skriver flera **magellan**-händelser i en loggfil (se tabellen nedan). En ny loggfil skapas en gång i veckan och lagras under

Windows 10: C:\Users\Public\Documents\Tecan\LogFiles\SystemAuditTrail

En kontrollsumma, som beräknas över hela loggfilen, används för att garantera loggfilens överensstämmelse och för att skydda filerna mot avsiktliga och oavsiktliga ändringar. Loggfilens överensstämmelse kontrolleras alltid vid start. Ogiltiga filer stängs omedelbart, markeras som ogiltiga genom att ".invalid" läggs till loggfilens namn, och en ny fil skapas i stället.

Gamla eller ogiltiga loggfiler flyttas till arkivmappen, där de skyddas mot radering.



## magellan-händelser som ska spåras

Händelse	Händelsetyp	Beskrivning
magellan Start	StartUp	magellan påbörjades.
magellan avslutas	Exit	magellan stängdes.
Lyckad inloggning av användare	Login	En användare har lyckats loggat in.
Misslyckad inloggning av användare	FailedLogin	Inloggningen misslyckades på grund av fel användarnamn eller lösenord.
Användaren loggar ut	Logout	En användare har loggat ut.
Ny arbetsyta	New	En ny arbetsyta skapades.
Ladda arbetsyta	Load	En befintlig arbetsyta laddades in i magellan.
Spara arbetsyta	Save	En arbetsyta har sparats.
Ny metod	New	En ny metod skapades.
Ladda metod	Load	En befintlig metodfil laddades in i magellan.
Spara metod	Save	En metod har sparats.
Ny prov-ID-lista	New	En ny prov-ID-lista har skapats.
Ladda prov-ID-lista	Load	En befintlig prov-ID-lista har laddats.
Spara prov-ID-lista	Save	En prov-ID-lista sparades.
Starta mätning	Measure	En ny mätning påbörjades.
Bifoga underskrift	Underskrift	En arbetsyta eller metodfil har signerats.
Exportera resultat	ExportASCII ExportExcel ExportASTM ExportSampleTra cking	Resultaten exporterades.
Resultat av utskrifter	Print	Resultaten skrevs ut.
Automatisk Arkivering	ArchiveFailed	Arkiveringen av en arbetsyta, metod eller prov-ID-lista misslyckades.
CRC-fel	CRCFailed	Kontrollsumman för den föregående loggfilen var ogiltig.
# 11. Beräkningar

# 11.1 Utvärdera resultaten – Beräkningsprocedur

Beräkningen följer nedanstående procedur beroende på inställningarna:

- 1. Precalculation (Förhandsberäkning)
  - 1.1 Spectra calculation (Spektraberäkning)
- 2. Check if data and evaluation settings fit together (Kontrollera om data och utvärderingsinställningar passar ihop)
- 3. Raw data statistic calculation (Beräkning av statistiska rådata)
- 4. Transformation calculation (Beräkning av transformation)
- 5. Kinetic parameter calculation (Beräkning av kinetiska parametrar)
- 6. Kinetic transformation calculation (Beräkning av kinetisk transformation)
- 7. Concentration calculation (Beräkning av koncentrationen)
- 8. Concentration transformation data (Koncentrationstransformationer)
- 9. Cutoff range determination (Fastställande av cutoff-intervall)
- 10.QC Validation (QC-Validering)
- 11. Statistic calculation (Beräkning av statistik)

Proceduren upprepas om

- QC-validering eliminerade värden med hjälp av elimineringsfunktionerna.
- Interpolationsparametrar i transformationerna krävs.

Lika värden för transformationer i kinetiska mätningar och mätningar med multipla inmärkningar kommer att undertryckas.

Vid beräkningar med olika spädningar kommer medelkoncentrationerna att avvisas.

Om utvärderingen avbryts på grund av fel kan de resultat som beräknats fram till dess visas.

Om vissa värden visar sig vara FALSE, ska följande dialogruta **Occurred Errors** (Fel inträffade) visas:

Textfältet <b>Message</b> (Meddelande)	Felmeddelandet och en kort förklaring visas i ett textfält.
Filknappon Sava as	Klicka nå knappan <b>Savo as filo</b> (Spara som fil) så kan

Filknappen Save as<br/>(Spara som)Klicka på knappen Save as file (Spara som fil) så kan<br/>felprotokollet sparas i en ASCII-fil.

Om kurvanpassningen misslyckades, beroende på användarens rättighet, visas en fråga där användaren kan besluta att ändra inställningarna för analystyp. Efter ändringarna börjar utvärderingsproceduren från början.

Om valideringen misslyckades visas, beroende på användarens rättighet, en ruta med ett felmeddelande eller fråga, där användaren kan besluta om han/hon vill fortsätta beräkningen ändå. Om beräkningen avbryts, gåt det inte att få tillgång cutoff-resultaten. På en utskrift kommer dock ett meddelande om misslyckad validering att finnas med.



Om värden eller inställningar har ändrats visas en meddelanderuta.

Message (meddelande)	Instrumentvärdena har ändrats. Vill du starta om beräkningen av resultatet?
Knappen <b>Yes</b> (Ja)	Klicka på knappen <b>Yes</b> (Ja) för att räkna om utvärderingsvärdena.
Knappen <b>No</b> (Nej)	Klicka på <b>No</b> (Nej) för att stänga meddelandet. Ingen ny beräkning kommer att göras.

# 11.1.1 Statistik

Om ett statistiskt värde beräknas över t.ex. n replikat med värden  $x_1...x_n$  av ett prov används följande formler:

Namn	Formel
Medelvärde, genomsnittligt värde	$\overline{x} = \frac{1}{n} \sum_{i=1}^{n} x_i$
Standardavvikelse	$s = \sqrt{\frac{1}{n-1} \sum_{i=1}^{n} (x_i - \bar{x})^2}$ och s=0 för n=1
Variationskoefficient i procent	$v = 100 \frac{s}{\overline{x}}$

Mer information finns i kapitel 11.3.5 Statistiska funktioner.

# 11.2 Reducering av spektradata

Spektradata innehåller mycket mer information än mätningar av enkla våglängder. Typisk information som extraheras från spektra är intensiteter vid specificerade våglängder eller vid hittade toppar. Under beräkningen är det också möjligt att beräkna spektrablankreducering och kvotberäkning.

# 11.2.1 Matematisk beskrivning

**Blankreducering** utförs genom att subtrahera blankspektrumet från provspektrumet.

**Utjämning** görs med Savitzky-Golay-algoritmen. Filterkoefficienterna beräknas med hjälp av ett polynom av grad 2. Utjämningsfaktorn behandlas som antalet punkter:

$$g_{i} = \sum_{n=-n_{L}}^{n_{R}} c_{n} f_{i+n}$$
där n<sub>R</sub>-n<sub>L</sub>+1=antal punkter

Utjämningsfönstret är symmetriskt, t.ex. för en utjämningsfaktor på 5 är nL=-2, nR=2.

För den nedre och övre gränsen av spektrumet är fönstret asymmetriskt, vilket innebär att  $n_L=0$ ,  $n_R=4$  eller  $n_L=-4$ ,  $n_R=0$ .

Före utjämning avlägsnas alla överflödsvärden från spektrumet.

#### Val av våglängd

• Pick wavelength (Välj våglängd) – Hämtar intensiteten för den valda våglängden



- Calculate ratio (Beräkna kvot) Beräknar kvoten genom att dividera intensiteterna vid de valda våglängderna
- Area Beräknar arean under spektrumet inom det valda våglängdsområdet
- Custom (Anpassad) Minskar spektrumet med hjälp av en användardefinierad formel

#### Normalize (Normalisera)

Reducerar spektrumet till ett intervall från 0 till 1.

#### Peak find (Hitta topp)

Spektrumet differentieras och alla nollpunkter i resultatet behandlas som potentiella toppar. Motsvarande toppar sorteras efter värde och alla ogiltiga toppar tas bort (t.ex. överskridanden, lokala minima, värden under det definierade tröskelvärdet). Det högsta värdet av de återstående elementeten väljs som toppvärde.

I praktiken är detta ofta det maximala värdet i spektrumet.

#### **Custom formulas (Anpassade formler)**

Möjlighet att definiera olika datareduktioner för spektra (se kapitel 11.3.8 Spektrafunctioner).



# VARNING ENSTAKA AVVIKELSER SOM ORSAKAS AV BRUS KOMMER OCKSÅ ATT UPPTÄCKAS SOM TOPPAR! DÄRFÖR ÄR DET STARKT REKOMMENDERAT ATT JÄMNA UT SPEKTRUMET INNAN DU ANVÄNDER PEAK FIND!

# 11.3 Hur man skriver en formel

# 11.3.1 Inledning

En mängd olika formler är nödvändiga för beräkningarna, cutoffs och valideringen. Dessa formler liknar syntaxen för BASIC-kommandon.

# Alla värden beräknas med dubbel precision även om de visas beroende på önskat talformat.

# 11.3.2 Formelvariabler

Alla tidigare definierade förkortningar för de identifierare som finns på analysplattan kan användas som variabler i beräkningar.

Till exempel:

PC1, SM1\_1, BL1...Medelvärden

PC1\_1, SM1\_1\_1\_1, BL1\_1...Individuella värden, individuella replikat



Anmärkning Programmet skiljer mellan små och stora bokstäver när det gäller dessa etiketter.

Det är också möjligt att arbeta med brunnskoordinater.

Till exempel: A2, B3, H12







i

Anmärkning Brunnskoordinater ska alltid skrivas med stora bokstäver.

Anmärkning Symbolen x hänvisar till det aktuella värdet i en brunn. Till exempel: x-BL1...subtraherar det tomma värdet (blanka) från det aktuella värdet.

Anmärkning De olika cyklerna för kinetiska mätningar kan kommas åt med index [ ], där "0" anger den första cykeln.

#### Till exempel:

SM1\_1[0]... Medelvärde för prov 1 i den första cykeln x[1].... värdet av den aktuella brunnen i den andra cykeln x[i] ... värdet av den aktuella brunnen och den aktuella cykeln

# Seriell kinetisk subtraktion – beräkning av skillnaden mellan kinetikcykler

Beräkningar av kinetiska data kan definieras genom att använda iterationsvariabeln "i" i formlerna. En numerisk härledning av en kinetisk kurva kan till exempel utföras med formeln x[i]-x[i-1], där varje kinetikcykel subtraheras med den föregående. Men om man använder denna formel skulle det automatiskt leda till ett fel för cykel 1. Ange därför formeln if(i>0) then(x[i]-x[i-1]) else ignore() i redigeringsfältet för transformation för att förhindra att den första cykeln subtraheras.

Användningen av konstanter i formler är identisk med alla andra variabler, den enda skillnaden är att en konstant kan innehålla tecken (t.ex. alfa, spädning).

Det finns två fördefinierade åtkomstmetoder.

# concX ...

återger koncentrationen av standarden i den aktuella brunnen

# dilX ...

återger spädningen av provet eller kontrollen av den aktuella brunnen

# Anmärkning

Om mer än ett set indata ska användas för beräkningar, välj lämpliga data från rullgardinsmenyn tillgängliga data, som aktiveras i redigeringsfälten för transformation på fliken Create/edit a method (Skapa/redigera en metod) (se kapitel 4.3.5 Transformed data (Transformerad data): Lägg till ny transformation och kapitel 7.4.9 Kontrollfält Transformerad data).



# Anmärkning

Genom att använda en av de matematiska funktioner som beskrivs i följande kapitel aktiveras en "intellisense" som stödjer korrekt skrivning av speciella formler.

Operator	Beskrivning
+	Plus
-	Minus
*	Multiplikation
1	Division
۸	Upphöjt till exponenten
	Symbolen som används för att representera uttrycket "upphöjt till" är ´^´. Detta tecken kan erhållas genom att trycka på SHIFT- tangenten tillsammans med tangenten nummer 6 (men inte på den numeriska knappsatsen)
	Brunnens värde kommer att ökas upphöjt till tre.
<	Mindre än
<=	Mindre än eller lika med
>	Större än
>=	Större än eller lika med
==	Lika med
!=	Är inte lika med
=	Tilldelad till

# Formeloperatorer



Anmärkning Skillnaden mellan lika (==) och tilldela till (=). Till exempel: Om (x == 0.000), så är x = 1.000 (Om det nuvarande värdet på x är lika med 0, sätt värdet på x till 1)



# 11.3.3 Formulelfunktioner

Uttrycken "och" och "eller" kan ingå i samma ekvation med hjälp av logiska funktioner. Resultatet av en logisk ekvation är antingen TRUE eller FALSE och är därför ett booleskt värde. De individuella ekvationerna måste omges av parenteser.

# Logiska uttryck

#### och

I ekvationer av detta slag blir resultatet TRUE om båda uttrycken stämmer, annars blir resultatet FALSE. Exempel:

Där värdet i brunnen är 0.3

(x>0.0) och (x<1) = TRUE

(x>0.0) och (x<0.1) = FALSE

# eller

I ekvationer av detta slag måste minst ett av de ingående uttrycken gälla för att man ska kunna få resultatet TRUE.

*Exempel:* där värdet i brunnen är 0.3

(x>0.0) eller (x<0.1) = TRUE (x>0.4) eller (x<0.1) = FALSK

# if(...) then(...) else(...)

Följande uttryck används i programmet och kan som exempel skrivas på följande sätt:

if (påstående) then Term A else Term B

Påstående:

Påståendet måste vara antingen sant eller falskt och skrivas inom parentes. Om påståendet görs med hjälp av logiska operatorer (och/eller) ska hela uttrycket skrivas här och omslutas av parenteser.

Termerna A och B:

Om påståendet är sant gäller term A och om påståendet är falskt gäller term B.

Det är inte alltid nödvändigt att inkludera uttrycket else. Om den inte finns med i ekvationen kommer inget svar att ges om ett felaktigt resultat uppstår.

Uttrycket "if(...) then(...) else(...)" kan också användas med booleska termer för QC-validering.

#### **Exempel:**

Kontrollera om värdet för en brunn ligger inom ett definierat intervall. Om resultatet är ja kommer brunnen att representeras av värdet 0.0, om resultatet är nej kommer värdet 1 att användas.

if (( x>-0.005) and (x<0.0))) then(0.0) else(1.0)



# 11.3.4 Grundläggande funktioner

# abs(argument)

	Denna funktion ger den absoluta summan av argumentet.
	Därför är (−1*x) där x<0 och (x) där x>= 0.
	Exempel:
	abs(-1) = 1
	abs(1) = 1
exp(argument)	
	Denna funktion gör att Eulervärdet (e) blir upphöjt till argumentet.
	<b>Exempel:</b> $exp(1) = e^1 = 2.718$
frac(argument)	
	Denna funktion separerar fraktionsdelen av argumentet.
	frac(): levererar bråkdelen av ett decimalvärde.
	<b>Exempel:</b> frac(1.7) = 0.7
int(argument)	
	Denna funktion separerar heltalsdelarna av argumentet.
	int(): levererar heltalsdelen av ett decimalvärde.
	<b>Exempel:</b> <i>int</i> (1.7) = 1
Log(argument)	
	log(arg1;arg2)
	Denna funktion tar logaritmen av arg2 till basen av arg1.
	Exempel:
	Brunnen har ett värde på 100.
	log(x;10)
	Här blir resultatet 0.5.
In(argument)	
	Denna funktion tar den naturliga logaritmen av argumentet.
	<b>Exempel:</b> <i>In</i> (10) = 2.303
lg(argument)	
	Denna funktion tar logaritmen med bas 10 av argumentet.
	<b>Exempel:</b> $lq(10) = 1$
round(argument)	
round(argument)	Denna funktion avrundar argumentet till ett heltalsvärde
	Exempel:
	För att avrunda 12 579 till två decimaler använder du följande formel:
	round(12.579*100)/100
	Resultatet är 12.58.

# sqr(argument)

Denna funktion gör att argumentet blir upphöjt till 2

**Exempel:** sqr(3) = 9

# sqrt(argument)

Denna funktion tar kvadratroten av argumentet

**Exempel:** sqrt(9) = 3

# 11.3.5 Statistiska funktioner



Anmärkning Använd enstaka data och inte genomsnitt eller medelvärde av data som indata för transformationer med statistiska funktioner.

# avg(argument) och mean(argument)

Denna funktion beräknar medelvärdet av ett argument. Argumentet måste vara en identifierare.

#### Exempel:

avg(SM1)

Medelvärdet av alla prover i den första försöksgruppen kommer att skapas. Dessutom kan medelvärdet av replikaten av ett värde beräknas:

avg(ST1\_1)

#### Medelvärden för alla replikat av den första standarden.

# median(argument)

Denna funktion bestämmer medianen för argumentet. Argumentet måste vara en identifierare. De individuella replikaten ordnas i storleksordning och medianen bestäms genom att ta det mittersta värdet i ordningen. Om antalet värden är ett jämnt antal bestäms medelvärdet av de två mittersta värdena.

# Exempel:

Median (NC1) Anta att NC1\_1=0.1, NC1\_2=0.05, NC1\_3=0.04 Medianvärdet för dessa negativa kontroller bestäms på följande sätt. Värdenas ordningsföljd ger: 0.04 0.05 0.1 Medianen är därför 0.05.

# medianPlate()

Medianen för hela plattan kan bestämmas med hjälp av denna funktion. De individuella värdena ordnas i storleksordning och medianen bestäms genom att ta det mittersta värdet i ordningen. Om antalet värden är ett jämnt antal bestäms medelvärdet av de två mittersta värdena.

#### Exempel:

medianPlate()

Anta att BL1=1, NC1=2, PC1=3, SM1\_1=4, SM1\_2=5 och inga andra brunnar definieras.

Medianen för hela plattan är 3.



# max(argument) och min(argument)

Denna funktion beräknar minimum/maximum för ett argument. Argumentet måste vara en identifierare.

#### Exempel:

max(SM1)

Maximala värdet av alla prover i den första försöksgruppen kommer att bildas.

min(ST1\_1)

Bestämmer miniminivån för de individuella replikaten av den första standarden.

#### maxAvg(argument) och minAvg(argument)

Denna funktion ger det maximala/minsta medelvärdet för argumentet. Argumentet måste vara en identifierare.

#### Exempel:

minAvg(SM1)

Om ett antal resultat hittas för SM1, kommer endast det lägsta genomsnittet att anges.

#### PointwiseCV(argument)

Den genomsnittliga punktvisa variationskoefficienten (CV) bestäms. Resultatet kan användas för att validera standardkurvan. Argumentet måste vara en identifierare.

#### Exempel:

PointwiseCV(ST1)

Anta

ST1\_1\_1=0.54 ST1\_1\_2=0.52 cv=2.668

ST1\_2\_1=0.72 ST1\_2\_2=0.77cv=4.746

ST1\_3\_1=1.08 ST1\_3\_2=0.99cv=6.148

Resultatet av pointwiseCV är genomsnittet av CV för ST1\_1, ST1\_2 och ST1\_3 och är 4.5209.

# Sum(argument)

Denna funktion beräknar summan av argumenten.

#### Exempel:

sum(SM1)

Standardavvikelsen för den negativa kontrollen i den första försöksgruppen kommer att bildas.

#### sum(ST1\_1)

Bestämmer summan av replikaten för den första standarden.

#### stddev(argument)

Denna funktion beräknar standardavvikelsen för ett argument. Argumentet måste vara en identifierare.

#### Exempel:

stddev(NC1)

Standardavvikelsen för den negativa kontrollen i den första försöksgruppen kommer att bildas.



# 11.3.6 Eliminationsfunktioner

Elimineringsfunktionerna kommer att användas vid valideringen. På så sätt kan användaren försäkra sig om att mätvärdena ligger inom ett giltigt intervall och radera eventuella felaktiga värden.

Tre olika elimineringsfunktioner kan användas när man definierar giltighetsintervallen. Den identifierare som ska påverkas av funktionen är alltid den första parametern som krävs när man definierar elimineringskriterierna.

Alla tre elimineringsfunktioner ger ett logiskt resultat. TRUE ges när det finns tillräckligt många giltiga mätvärden tillgängliga och FALSE ges när färre än vad som krävs är giltiga.

Ogiltiga mätvärden markeras med ett utropstecken.

# eliminate (arg1; arg2; arg3; arg4)

Valideringsintervallet definieras med hjälp av ett definierat värde.

Denna funktion eliminerar värden som ligger utanför det valideringsområde som definieras i arg2.

arg1: Identifieringsnamn (NC1,PC1).

arg2: Värde som används för att representera intervallet, beroende på arg4. arg3: Antal giltiga, individuella värden som måste vara tillgängliga för att ett medelvärde ska kunna genereras. Om för få värden finns tillgängliga efter elimineringen får du resultatet FALSE.



Anmärkning För arg3: ett värde på "2" eller högre måste definieras. "1" accepteras inte.

arg4: Representerar valet om testningen ska använda intervallet averagearg2 till average+arg2 (argument 1) eller om det ska kontrolleras om de individuella värdena ligger under average+arg2 (argument 0). I detta fall (argumentet 0) finns det ingen nedre gräns.

Elimineringsfunktionen bestämmer här genomsnittet för arg1. Det värde som har den största absoluta avvikelsen från medelvärdet kontrolleras sedan för att säkerställa att det ligger inom intervallet enligt arg4. Om värdet ligger utanför detta intervall betraktas det som ogiltigt. I detta fall börjar genomsnittsberäkningen på nytt, utan att ta hänsyn till det tidigare upptäckta

ogiltiga värdet, och fortsätter på detta sätt tills inga fler felaktiga värden kan hittas.

Resultatet TRUE kommer att levereras när tillräckligt många giltiga individuella värden finns tillgängliga.

# Exempel:

eliminate(NC1;0.15;2;0)

Medelvärdet kommer att genereras för de negativa kontrollerna. Därefter kan man se om de individuella negativa kontrollerna ligger över NC1+0.15. (Det finns ingen nedre gräns). Om så är fallet ska det/de felaktiga värdet/värdena elimineras på det sätt som beskrivs ovan. Om minst två individuella värden är giltiga vid slutförandet returneras resultatet TRUE, annars returneras resultatet FALSE.

# eliminatePerc (arg1; arg2; arg3; arg4)

Valideringsintervallet definieras med hjälp av en procentsats av medelvärdet.

Denna funktion eliminerar värden på samma sätt som beskrivs för funktionen eliminera. Valideringsområdet definieras med hjälp av en procentandel (arg2) av identifieraren.



# Exempel:

elimineraPerc(NC1;10;2;1)

Individuella värden är ogiltiga om de avviker med mer än 10 % under eller över genomsnittet. Vid slutförandet ska minst två individuella värden vara giltiga för att de negativa kontrollerna skall förklaras vara giltiga.

# eliminateRange (arg1; arg2; arg3; arg4)

Denna funktion eliminerar värden som ligger utanför det valideringsområde som definieras av argumenten 2 och 3.

arg1: Identifieringsnamn (NC1,PC1).

arg2: Lägre gräns för det valda intervallet

arg3: Övre gräns för det valda intervallet

arg4: Antalet giltiga, individuella värden som måste finnas för att fortsätta. Om för få värden finns tillgängliga efter elimineringen får du resultatet FALSE.

Resultatet TRUE kommer att levereras när tillräckligt många giltiga individuella värden finns tillgängliga.

# Exempel:

eliminateRange(NC1; 0.0; 0.1; 2)

De individuella värdena för de negativa kontrollerna måste ligga inom intervallet 0.0 och 0.1 för att vara giltiga. För att uppnå resultatet TRUE måste minst två av värdena anses vara giltiga.

# eliminateCV (arg1; arg2; arg3)

Denna funktion eliminerar replikaten tills CV för de återstående replikaten är lägre än det givna CV (arg2). Replikaten elimineras stegvis, med början med det värde som har den största skillnaden mot medelvärdet.

arg1: Identifieringsnamn

arg2: Gränsen för det accepterade CV-värdet

arg3: Antalet giltiga, individuella värden som måste lämnas kvar för att ge ett positivt resultat. Om för få värden finns tillgängliga efter elimineringen får du resultatet FALSE.

Resultatet TRUE kommer att levereras när tillräckligt många giltiga individuella värden finns tillgängliga.

#### Exempel:

eliminateCV(NC1; 15; 3)

Det beräknade CV-talet för de individuella värdena för den negativa kontrollen ska vara lägre än 15 %. Efter elimineringen måste det finnas minst 3 replikat kvar för att få TRUE som resultat.

# countDeleted(arg1; arg2)

Den här funktionen kontrollerar om det finns tillräckligt många giltiga värden och returnerar TRUE eller FALSE.

arg1: Den identifierare vars kvantitet ska kontrolleras.

arg2: Det lägsta antalet replikat som måste vara tillgängligt.

#### Exempel:

countDeleted(NC1; 2)

Den här funktionen fungerar tillsammans med funktionerna Eliminera i ekvationer:

Valideringen innehåller följande rad:



om (NC1>0.5) elimineraRange(NC1; 0; 0.5; 2

Om medelvärdet för den negativa kontrollen ligger över 0.5 ska alla replikat som ligger utanför intervallet utelämnas.

Denna eliminering kommer att behandlas en gång. Om värden för eliminering hittas, görs en ny beräkning för att se till att genomsnittet ligger under 0.5 och resultatet blir TRUE.

I detta skede är det ännu inte säkert att det nödvändiga antalet replikat finns tillgängliga. Detta bestäms sedan med hjälp av countDeleted-funktionen i ett elsemeddelande:

om (NC1>0.5) elimineraRange(NC1; 0; 0.5; 2) annars countDeleted(NC1; 2)

I ekvationsprocessen (i den andra körningen efter elimineringen) används elseangivelsen. Funktionen countDeleted kontrollerar sedan om den givna identifieraren i arg1 finns tillgänglig i den mängd som anges i arg2. Om detta är fallet får du resultatet TRUE, annars FALSE.

#### Steg för steg exempel eliminera respektive elimineraPerc

Den enda skillnaden mellan eliminate och eliminatePerc är sättet att beräkna gränserna för det giltiga området (se ovan). Beräkningssekvensen är densamma för båda funktionerna. För den stegvisa beskrivningen används eliminering.

#### eliminate(NC1;0.15;2;0)

Steg 1: De individuella värdena sorteras efter sitt avstånd till genomsnittet.

$ NC1_1 - \overline{NC1} $	0.217-0.288	0.071		NC1_2 = 0.439
$ NC1_2 - \overline{NC1} $	0.439-0.288	0.151	→	NC1_3 = 0.208
$ NC1_3 - \overline{NC1} $	0.208-0.288	0.08		NC1_1 (0.217)

Steg 2: Beräkning av den övre gränsen: 0.288+0.15 = 0.438

Steg 3: första jämförelsen:

0.439 > 0.438 och därmed utanför det giltiga intervallet  $\rightarrow$  Steg 4: Värdet för NC1\_2 elimineras och medelvärdet och gränsen beräknas på nytt.

Steg 4: Nytt genomsnitt = (0.217+0.208)/2 = 0.2125

Ny gräns = 0.2125+0.15 = 0.3625

Steg 5: andra jämförelsen:

0.208 <= 0.3625 → giltigt

0.217 <= 0.3625 → giltigt

#### Resultat:

Replikatet NC1\_2 passar inte in i det givna intervallet och har därför eliminerats.

Funktionen returnerar TRUE eftersom det fortfarande finns två replikat kvar.

Anmärkning: I varje beräkningscykel (steg 1 till steg 4) raderas ett replikat bort (det replikat som har det största avståndet till genomsnittet). Cykeln upprepas tills inget replikat har raderats under jämförelsesteget i cykeln.



# 11.3.7 Andra funktioner

# ignore()

Denna funktion utelämnar värdena för valda brunnar så att de inte tas med i beräkningarna. Ignorerade värden visas med ett inledande '!' och hanteras som eliminerade värden.

# Exempel:

if(x<0.9\*SM1\_1) then ignore()

# isInvalid()

Den här funktionen ger ett booleskt värde som anger om värdena i brunnen är giltiga eller inte. Ogiltiga värden kan till exempel orsakas av att läsaren har för höga värden.

#### Exempel:

if(isInvalid()) då x=3.0

Om brunnen innehåller ett ogiltigt värde tilldelas värdet 3.0. På så sätt är det möjligt att göra ytterligare beräkningar trots sådana värden.

# calcAlways(argument)

Denna funktion gör det möjligt att beräkna formler på brunnar oberoende av värdet. Argumentet kan vara en giltig formel. Formeln beräknas även om värdet är dolt eller ogiltigt. Denna funktion kan användas för att beräkna resultat som inte beror på värdet av den använda brunnen.

# Exempel:

calcAlways(A)

Värdet för parameter A i formeln för standardkurvan returneras.

concX

Denna funktion kan användas vid koncentrationstransformationer. Resultatet av **concX** är brunnens ursprungliga koncentration och kan därför endast användas i brunnar med standarder som identifierare.

# Exempel:

x – concX anger skillnaden mellan den beräknade och den ursprungliga koncentrationen i brunnen.

# 11.3.8 Spektrafunctioner

# smooth(numPts)

Utjämnar med det angivna antalet punkter.

minimum()

Levererar det minsta värdet som ett datapar (våglängd, intensitet).

maximum()

Levererar det maximala som datapar (våglängd, intensitet).

# normalize(wl)

Normaliserar genom att använda intensiteten vid den givna våglängden.



# peak(threshold;wlStart;wlEnd;peakIndex:percLimit)

Levererar toppen som ett dataset (intensitet, våglängd, bredd, area). Toppar som är lägre än tröskelvärdet kommer inte att hittas.

Sökning av toppar görs i det givna våglängdsområdet.

peakIndex = -1  $\rightarrow$  återger max. topp

peakIndex =  $-2 \rightarrow$ återger min. topp

peakIndex = 0, 1, 2, ...  $\rightarrow$  returnerar toppen vid det angivna indexet

percLimit är en procentsats som används för att beräkna bredd och area.

# numPeaks(threshold;wlStart;wlEnd)

Ger antalet toppar som hittats i det givna våglängdsområdet. Toppar som är lägre än tröskelvärdet kommer inte att hittas.

# derive(degree)

Beräknar derivatet av den givna graden. Giltiga grader är 1 och 2, t.ex. derive(2).

# intensity(function)

Ger intensiteten för antingen ett datapar eller en topp; t.ex. intensity(minimum()).

# wavelength(function)

Levererar våglängden för antingen ett datapar eller en topp, t.ex. våglängd(maximum())

# width(function)

Ger bredden av en topp, t.ex. width(peak0.5;450;650;-1;50)).

# area(function)

Ger område av en topp, t.ex. area(peak0.5;450;650;-1;50)).

# 11.3.9 Exempel

# Transformationer

Transformationer ändrar brunnens aktuella värde. Resultatet måste alltid vara ett numeriskt värde.

#### Blank reduction (Blankreducering):

x-BL1

Det tomma värdet (blank) kommer att dras av från brunnens nuvarande värde.

#### Blanking of a kinetic run (Blankreducering av kinetisk omgång:

x-x[0]

Värdet för den första punkten i kinetiken används som blank och kommer att härledas från värdena för alla kinetikcykler.



Anmärkning Indexet "0" representerar den första mätningen av en kinetisk omgång.



# Ratio (Kvot):

x/PC1 \* 100

Beräknar kvoten mellan ett prov och en kontroll i procent.

#### **DNA/Protein ratio:**

'Label1'!x/'Label2'!x

För att beräkna kvoten mellan DNA och protein måste en mätning med multipla inmärkningar definieras. Vid den första mätningen används en våglängd på 260 nm och vid den andra mätningen en våglängd på 280 nm.

#### Conditional result (Villkorligt resultat)

Om (x<0.0) då 0 annars x

Om värdet av brunnen är mindre än 0 representeras det av 0, annars förblir det oförändrat.

# Cutoff

Cutoff anger vilka gränser som ska tillämpas vid utvärderingen. Resultatet måste alltid vara ett numeriskt värde.

# Märkning: POS

Gräns: NC1\*1.15

Märkning: ??

Gräns: NC1\*0.85

# Märkning: -

Resultatet är positivt (POS) när absorbansen är mer än 15 % högre än genomsnittet av de negativa kontrollerna och negativt (-) när den är mer än 15 % lägre än genomsnittet. Ett resultat anges som öppet (???) om det ligger mellan dessa två värden.

# **QC Validation (QC-validering)**

Valideringen syftar till att undersöka testets giltighet. Resultatet måste alltid vara ett logiskt uttalande, TRUE eller FALSE.

# (BL1>-0.005) och (BL1<0.120)

Om medelvärdet av det tomma värdet ligger inom intervallet -0.005 till 0.12, kommer resultatet att vara TRUE.

# eliminateRange(NC1;-0.005;0.12;2)

Om minst två negativa kontroller ligger inom intervallet -0.005 till 0.12, kommer resultatet att vara TRUE.

# abs(PC1-NC1)>0.25

Avståndet mellan medelvärdena för de positiva kontrollerna och de negativa kontrollerna skall vara större än 0.25. Om det är så kommer resultatet att vara TRUE.

# **11.4** Typer av analys av standardkurvor

# 11.4.1 Definitioner

Givet n baspunkter  $(x_1, y_1), ..., (x_n, y_n), x_1 < ... < x_n$ 

Namn	Formel
Correlation coefficient (korrelationskoefficient)	$r = \frac{\sum_{i=1}^{n} (x_i - \bar{x})(y_i - \bar{y})}{\sqrt{\sum_{i=1}^{n} (x_i - \bar{x})^2 \sum_{i=1}^{n} (y_i - \bar{y})^2}}$
Average square deviation (genomsnittlig kvadratavvikelse)	$d = \sqrt{\frac{1}{n} \sum_{i=1}^{n} (y_i - f(x_i))^2}$
Goodness of fit (anpassningsgrad)	goodness = 1 - d

# 11.4.2 Parametrar för analystyp

Parametern för analystypen (t.ex. lutningen A och skärningspunkten B för linjär regression) kan användas i beräkningen. De tillgängliga parametrarna beskrivs nedan. **Om det finns mer än en försöksgrupp** har parametrarna för analystypen suffixet "B" för den andra försöksgruppen, "C" för den tredje försöksgruppen och så vidare. Exempel: Lutningen för en linjär regression i försöksgrupp 3 kan anges som "AC".

# 11.4.3 Felmeddelanden

Om något av värdena i de beräknade koncentrationerna ligger utanför det intervall som anges i standarderna, ska följande inmatning **>Max** eller **<Min** visas i plattans brunnar. Om användaren ändå vill ta med dessa värden i beräkningen måste alternativet Extrapolation (Extrapolering) väljas.

Om kurvan inte är strikt monoton genereras ett felmeddelande.

Om det finns flera lösningar för en koncentrationsberäkning för ett uppmätt värde visas inmatningen MultPt i plattans brunnar.

# 11.4.4 Point to point (Punkt-till-punkt)

Med den här processen sammanfogas de intilliggande baspunkterna med hjälp av en rät linje.

Minst 2 baspunkter krävs för denna beräkning.

Extrapolering är inte möjlig. Det finns inga parametrar som kan användas som variabler i transformationerna.

Om ingen koncentration kan beräknas visas NoCalc i plattans brunnar.

#### Matematisk beskrivning

Interpoleringsfunktion:

$$f:[x_1, x_n] \to R$$
  
$$x \mapsto y_i + \frac{y_{i+1} - y_i}{x_{i+1} - x_i} \cdot (x - x_i) \quad \text{if } x \in [x_i, x_{i+1}]$$



# 11.4.5 Linear Regression (Linjär regression)

En rak linje bildas där summan av de kvadratiska avvikelserna för alla baspunkter är ett minimum.

Minst 2 baspunkter krävs för denna beräkning.

Vid behov kan extrapolering också genomföras i denna process. I en utvärdering med en försöksgrupp fungerar lutningen och y-axelns skärningspunkt som variablerna A och B i transformationerna. Korrelationskoefficienten r kan också användas.

I dialogrutan Standardkurva visas följande ytterligare information:

Formel med koefficientvärdena A (lutning) och B (skärningspunkt)

Genomsnittlig kvadratisk avvikelse d

Absoluta värdet av korrelationskoefficienten r

#### Matematisk beskrivning

Approximationsfunktion:

$$f:[x_1,x_n]\to R$$

$$x \mapsto A \cdot x + B$$

där A och B bestäms genom att minimera felfunktionen

$$err(A, B) = \sum_{i=1}^{n} (f(x_i) - y_i)^2$$

Lösningen är unik om

$$rank \begin{pmatrix} x_1 & 1 \\ \vdots & \vdots \\ x_n & 1 \end{pmatrix} = n$$
, vilket är sant om  $x_i \neq x_j \forall i, j = 1,...,n$  (se det allmänna

villkoret)

# 11.4.6 Non-Linear Regression (Icke-linjär regression)

Denna approximation är särskilt utformad för hyperboliska data. Idealt är de uppmätta värdena för låga koncentrationer höga och de uppmätta värdena för höga koncentrationer låga och det intresseområdet ligger i det höga koncentrationsområdet.

Den parameterberäkning som stöds av den linjära regressionen x/y används för y.

En linjär regression utförs med hjälp av de transformerade data.

Resultatet uttrycks som lutningen k och skärningspunkten d.

Utifrån detta resultat kan A = 1/k och B = d/k bestämmas.

Minst 2 baspunkter krävs för denna beräkning.

Vid behov kan extrapolering också genomföras i denna process. I en utvärdering med en försöksgrupp fungerar koefficienterna A och B som variabler i transformationerna. Korrelationskoefficienten r kan också användas. Approximationen tillämpas alltid på de oskalade baspunktsvärdena.

I dialogrutan Standardkurva visas följande ytterligare information:

Formel med koefficientvärden A och B

Genomsnittlig kvadratisk avvikelse d

Absoluta värdet av korrelationskoefficienten r





# VARNING

# DENNA APPROXIMATION ÄR INTE KONTINUERLIG OCH DÄRFÖR KAN FELAKTIGA DATA ORSAKA STORA AVVIKELSER MELLAN DE GIVNA BASPUNKTERNA OCH DE BERÄKNADE PUNKTERNA!

# Matematisk beskrivning

Approximationsfunktion:

$$f:[x_1, x_n] - \{-b\} \to R$$
$$x \mapsto \frac{A \cdot x}{B + x}$$

där A och B bestäms genom att lösa det linjära regressionsproblemet för de transformerade baspunkterna

$$(x_1, \frac{x_1}{y_1}), \dots, (x_n, \frac{x_n}{y_n})$$

Linjär regression:

$$g:[x_1,x_n] \to R$$

 $x \mapsto k \cdot x + d$  , som minimerar

$$err(k,d) = \sum_{i=1}^{n} (g(x_i) - \frac{x_i}{y_i})^2$$

Parametrarna A och B beräknas från k och d genom

$$A = \frac{1}{k} \quad B = \frac{d}{k}$$

Denna funktion f är inte kontinuerlig vid –B.

# 11.4.7 Polynomial (Polynom)

2:a grad: parabolisk eller kvadratisk

3:e grad: kubisk

Användaren kan välja mellan en beräkning av ett polynom av 2:a graden eller ett polynom av 3:e graden.

Minst 3 baspunkter krävs för beräkning av ett polynom av 2:a graden, medan minst 4 krävs för beräkning av ett polynom av 3:e graden.

Vid behov kan extrapolering också genomföras i denna process. Parametrarna A =  $a_0 B = a_1$  och C =  $a_2$  kan användas i transformationer med polynom av 2:a graden. För polynom av 3:e graden, kan variablerna A =  $a_0$ , B =  $a_1$ , C =  $a_2$  och D =  $a_3$  genomföras.

I dialogrutan Standardkurva visas följande ytterligare information:

Formel med koefficientvärdena A, B och C för polynom av 2:a graden eller formel med koefficientvärdena A, B, C och D för polynom av 2:a graden

Genomsnittlig kvadratisk avvikelse



#### Matematisk beskrivning

Approximationsfunktion:

$$\begin{split} &f:[x_1,x_n] \to R \\ &x \mapsto \sum_{i=0}^{order} a_i \cdot x^i \\ &\text{, grad= 2 eller 3} \end{split}$$

(n>grad)

där engrad,..., ao bestäms genom att minimera felfunktionen

$$err(a_{order},...,a_0) = \sum_{i=1}^n (f(x_i) - y_i)^2$$

Lösningen är unik om

$$rank \begin{pmatrix} x_1^{order} & \dots & x_1 & 1 \\ \vdots & \vdots & \vdots & \vdots \\ x_n^{order} & \dots & x_n & 1 \end{pmatrix} = order + 1$$
, vilket är sant om  $x_i \neq x_j \forall i, j = 1, \dots, n$ 

# 11.4.8 Kubisk spline

De intilliggande baspunkterna kommer att sammanfogas genom beräkning av ett polynom av 3:de graden. Parametrarna bestäms med hjälp av villkoret "not-a-knot".

Minst 3 baspunkter krävs för denna beräkning.

Extrapolering är inte möjlig. Det finns inga parametrar som kan användas som variabler i transformationerna.

#### Matematisk beskrivning

Funktioner för interpolering:

Styckvis polynom av 3:e graden.

Kontinuerlig andra härledning vid alla baspunkter.

$$\int_{x_1}^{x_n} \left(\frac{d^2 f}{dx^2}\right)^2 dx$$
 Minimerar integralen

, som är en förenklad term för

deformationsenergin för en spline.

Villkoret "not-a-knot": vid  $x_2$  och  $x_{n-1}$  är även den tredje härledningen kontinuerlig.

# 11.4.9 Akima

Akima-interpolationen skapar i vissa fall en mjukare kurva än splineinterpolationen.

Minst 3 baspunkter krävs för denna beräkning.

Extrapolering är inte möjlig. Det finns inga parametrar som kan användas som variabler i transformationerna.

#### Matematisk beskrivning

Denna interpoleringsmetod använder styckvisa polynomer av 3:e graden.

Det styckvisa interpolationspolynomet och dess första härledning är kontinuerliga.



# Method (Metod)

Lutningarna  $m_i$  i alla baspunkter uppskattas från grannpunkterna med hjälp av en särskild funktion:

$$q_i = \frac{y_i - y_{i-1}}{x - x}$$

 $x_i - x_{i-1}$  är lutningarna för den linjära interpolatorn mellan punkt i och punkt i-1, i=2,...,n

$$m_{i} = \frac{q_{i} |q_{i+2} - q_{i+1}| + q_{i+1} |q_{i} - q_{i-1}|}{|q_{i+2} - q_{i+1}| + |q_{i} - q_{i-1}|}, i=3,...,n-2$$

Särskilda fall:

 $\operatorname{Ar} \mathbf{q}_{i-1} = \mathbf{q}_i \operatorname{och} \mathbf{q}_{i+1} = \mathbf{q}_{i+2}$ , ange

Är  $\mathbf{q}_i = \mathbf{q}_{i+1}$ , ange  $\mathbf{p}'$  ( $\mathbf{x}_i$ ) =  $\mathbf{q}_i = \mathbf{q}_{i+1}$ . Är  $\mathbf{q}_{i-1} = \mathbf{q}_i$  och  $\mathbf{q}_{i+1} \Box \mathbf{q}_{i+2}$ , så är  $\mathbf{y}'_i = \mathbf{q}_i$  (analogt för  $\mathbf{q}_{i+1} = \mathbf{q}_{i+2}$ ).

$$m_i = \frac{q_i + q_{i+1}}{2}$$

För indexen 1, 2, n-1, n kan lutningarna inte uppskattas med denna algoritm.

Nu har vi följande villkor för den 3:e gradens interpolationspolynom pi, i=1, ..., n-1

pi'(xi) =mi

pi'(x<sub>i+1</sub>) =m<sub>i+1</sub>

pi'(xi) =mi

 $p_i'(x_{i+1}) = m_{i+1}$ 

vilket är fyra villkor för varje 3:e gradens interpolationspolynom pi.

# 11.4.10 LogitLog

LogitLog är utformad för sigmoidala standardkurvor. LogitLog ger alltid en Sformad kurva. Kurvan kommer asymptotiskt att nå en övre och en undre gräns, som oberoende av varandra kan fastställas som minsta och maximala gränser. Om användaren inte har fastställt några gränser här används det minsta och maximala baspunktsvärdena. Approximationen tillämpas alltid på de oskalade baspunktsvärdena.

Minst 4 baspunkter krävs för denna beräkning.

Vid behov kan extrapolering också genomföras i denna process. Parametrarna A, B, C och D kan användas som parametrar i transformationerna.

I dialogrutan Standardkurva visas följande ytterligare information:

Koefficientvärden A, B, C och D

Genomsnittlig kvadratisk avvikelse d

Korrelationskoefficientens absoluta värde r (korrelation mellan data och anpassning)

#### Restriktioner för processtillämpning

LogitLog och approximation med fyra parametrar får endast användas när man arbetar med standardkurvor som uppfyller följande krav:

- Sigmoid funktion tillämpad på standarddata i förhållande till koncentrationerna
- Symmetriskt genererad kurva i förhållande till IC50-värdet



 För att beräkna kurvan korrekt är det nödvändigt att ha ett koncentrationsvärde på noll och ett värde på oändlig standardkoncentration (mättat värde), så att asymptoterna kan bestämmas. Om dessa två värden inte finns i standardkurvan kan de definieras manuellt med hjälp av alternativet More standard curve parameters (Fler parametrar för standardkurvor) (Min, Max). Om inga Min- och Max-värden har definierats, kommer magellan basera processen på de minsta respektive största värdena.



VARNING OM DESSA KRAV INTE UPPFYLLS KAN APPROXIMATION AV DENNA TYP INTE ANVÄNDAS KORREKT.

# Matematisk beskrivning

LogitLog-regressionen är en funktion

$$f:[x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{1 + \left(\frac{x}{C}\right)^B}$$

för beskrivning av sigmoid korrelation av data.

Parametern kan tolkas som:

$$A = \lim_{x \to 0} f(x)$$
$$D = \lim_{x \to \infty} f(x)$$
$$f(C) = \frac{A + D}{2}$$

A, D bestäms som minimum respektive maximum (eller vice versa om funktionen är avtagande).

Därefter löses problemet med linjär regression för transformerade baspunkter.

$$X = \log_{10} x$$
$$Y = \ln \left( \frac{\frac{D - y}{D - A}}{1 - \frac{D - y}{D - A}} \right)$$

 $g: [X_i, Y_i] \to R$  $x \mapsto k \cdot x + d$  , som minimerar felfunktionen

$$err(k,d) = \sum_{i=1}^{n} (g(x_i) - Y_i)^2$$

Parametrarna B och C bestäms utifrån k, d:

$$B = -m \cdot \log_{10}(e)$$
$$C = e^{\frac{b}{B}}$$



Parametrarna A,B,C,D har tilltalande geometriska tolkningar.

A = respons vid en dos (x-värde) på noll (tänk på att x=0 inte visas i logaritmiska diagram)

D = svar vid en oändligt hög dos

C = är symmetripunkten på kurvan (IC50), och under denna punkt är kurvan en spegelbild

B = är ett mått på hur brant kurvan är vid inflexionspunkten; i själva verket är det (med en möjlig teckenändring) kurvans lutning i logit-form

Observera att kurvan alltid är avtagande (om A>D) eller ökande (om A<D).

# 11.4.11 Fyra parametrar

Approximationen med fem parametrar måste i princip uppfylla samma krav på baspunkter som LogitLog.

Den resulterande kurvan kommer att tas fram med hjälp av Nelder och Mead Downhill-Simplex-metoden. Denna metod ger en allt mer exakt approximation av resultatet under hela intervallet. Metoden med fyra parametrar kräver betydligt mer tid för att behandlas. Precis som med LogitLog skapas en S-formad kurva.

Minst fyra baspunkter krävs för denna beräkning.

Vid behov kan extrapolering också genomföras i denna process. Parametrarna A, B, C och D kan användas som parametrar i transformationerna. Approximationen tillämpas alltid på de oskalade baspunktsvärdena.

I dialogrutan Standardkurva visas följande ytterligare information:

Koefficientvärden A, B, C och D

Genomsnittlig kvadratisk avvikelse d

Korrelationskoefficientens absoluta värde r (korrelation mellan data och anpassning)

#### Restriktioner för processtillämpning

Se kapitel 11.4 Typer av analys av standardkurvor – LogitLog för ytterligare information.

#### Matematisk beskrivning

Approximationsfunktion:

$$f:[x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{1 + \left(\frac{x}{C}\right)^B}$$

#### Method (Metod)

Först beräknas en LogitLog-approximation.

Parametrarna A, B, C och D optimeras med hjälp av Downhill-Simplex-algoritmen, en minimeringsalgoritm som endast kräver funktionsutvärderingar men ingen funktionshärledning.

Algoritmen stannar om en noggrannhet på 0.001 uppnås (lyckad) eller om det maximala antalet 10 000 iterationer överskrids (misslyckande) innan den givna noggrannheten uppnås.

#### Bibliografi

Nelder, J.A., and Mead, R. 1965, Computer Journal, vol. 7, pp 308-313



# 11.4.12 Fyra parametrar – Marquardt

Approximationen med fem parametrar måste i princip uppfylla samma krav på baspunkter som LogitLog.

Den resulterande kurvan kommer att tas fram med hjälp av Levenberg-Marquardt-metoden. Denna metod ger en allt mer exakt approximation av resultatet under hela intervallet. Metoden med fyra parametrar kräver betydligt mer tid för att behandlas. Precis som med LogitLog skapas en S-formad kurva.

Minst fyra baspunkter krävs för denna beräkning.

Vid behov kan extrapolering också genomföras i denna process. Parametrarna A, B, C och D kan användas som parametrar i transformationerna. Approximationen tillämpas alltid på de oskalade baspunktsvärdena.

I dialogrutan Standardkurva visas följande ytterligare information:

- Koefficientvärden A, B, C och D
- Genomsnittlig kvadratisk avvikelse d
- Korrelationskoefficientens absoluta värde r (korrelation mellan data och anpassning)

#### Restriktioner för processtillämpning

Se kapitel 11.4 Typer av analys av standardkurvor – LogitLog för ytterligare information.

#### Matematisk beskrivning

Approximationsfunktion:

$$f:[x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{1 + \left(\frac{x}{C}\right)^B}$$

# Method (Metod)

Först beräknas en LogitLog-approximation.

Parametrarna A, B, C och D optimeras med Levenberg-Marquardt-algoritmen , en iterativ teknik som hittar ett lokalt minimum för en funktion som uttrycks som summan av kvadrater av icke-linjära funktioner.

Algoritmen stannar om en noggrannhet på 1E-7 (FLT\_EPSILON) uppnås (lyckad) eller om det maximala antalet 30 000 iterationer överskrids (misslyckande) innan den givna noggrannheten uppnås.

# 11.4.13 Fem parametrar – Marquardt

Approximationen med fem parametrar måste i princip uppfylla samma krav på baspunkter som LogitLog. Den sigmoidala kurvan kan dock vara osymmetrisk.

Den resulterande kurvan kommer att tas fram med hjälp av Levenberg-Marquardt-metoden Denna metod ger en allt mer exakt approximation av resultatet under hela intervallet. Metoden med fem parametrar kräver betydligt mer tid för att behandlas. Precis som med LogitLog skapas en S-formad kurva.

Minst fem baspunkter krävs för denna beräkning.

Vid behov kan extrapolering också genomföras i denna process. Parametrarna A, B, C, D och E kan användas som parametrar i transformationerna. Approximationen tillämpas alltid på de oskalade baspunktsvärdena.



I dialogrutan Standardkurva visas följande ytterligare information:

- Koefficientvärden A, B, C, D och E
- Genomsnittlig kvadratisk avvikelse d
- Korrelationskoefficientens absoluta värde r (korrelation mellan data och anpassning)

#### Restriktioner för processtillämpning

Se kapitel 11.4 Typer av analys av standardkurvor – LogitLog för ytterligare information.

#### Matematisk beskrivning

Approximationsfunktion:

$$f: [x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{(1 + (\frac{x}{C})^B)^E}$$

#### Method (Metod)

Parametrarna A, B, C, D och E optimeras med Levenberg-Marquardt-algoritmen, en iterativ teknik som hittar ett lokalt minimum för en funktion som uttrycks som summan av kvadrater av icke-linjära funktioner.

Algoritmen stoppas om en noggrannhet på 1E-7 (FLT\_EPSILON) uppnås (lyckad) eller om det maximala antalet 30 000 iterationer överskrids (misslyckad) innan den givna noggrannheten uppnås.

# 11.4.14 Viktning för fyra/fem parametrar – Marquardt/Polynomial Fit

Viktning påverkar standardoptimeringsalgoritmen.

Algoritmen försöker optimera den viktade summan av de minsta kvadraterna.

En viktningsfaktor på 1 innebär ingen viktning.

Viktningsfaktorer som är större än ett innebär att den angivna punkten har högre prioritet för den resulterande anpassningen.

Viktningsfaktorer som är mindre än ett anger att punkten har lägre prioritet.



# VARNING

VIKTER BÖR ENDAST ANVÄNDAS OM DET FINNS SERIÖSA STATISTISKA BEVIS FÖR ATT DE INSAMLADE DATA OCH ALGORITMEN ÄR LÄMPLIGA.



# Matematisk beskrivning

Genomsnittlig viktade kvadratisk avvikelse

$$d = \sqrt{\frac{1}{\sum_{i=1}^{n} k_i} \sum_{i=1}^{n} k_i (y_i - f(x_i))^2}$$

#### Automatisk beräkning av vikter med hjälp av varians:

Vikterna beräknas automatiskt med 1/SD<sup>2</sup> där SD är standardavvikelsen för replikaten av den faktiska baspunkten.

Detta innebär att data med hög standardavvikelse viktas mindre än data med låg standardavvikelse.



# VARNING VIKTNING MED HJÄLP AV VARIANS ÄR FARLIG NÄR FÖR FÅ REPLIKAT ANVÄNDS, EFTERSOM STANDARDAVVIKELSEN DÅ KAN HA EN HÖG SLUMPMÄSSIG ANDEL SOM AVSEVÄRT PÅVERKAR OPTIMERINGSALGORITMEN.

Automatisk beräkning av vikter med hjälp av relativa vikter:

Vikterna beräknas automatiskt med 1/Y<sup>2</sup> där Y är medelvärdet för den faktiska baspunkten. Detta innebär att algoritmen minimerar de relativa avstånden mellan data och kurvan.

# 11.5 Beräkning i spädningsserien

# 11.5.1 Upptäckt av spädningsserier

Följande kriterier måste vara uppfyllda för att tvinga fram **magellan** att upptäcka en spädningsserie på layouten:

- Prov med minst fyra replikat
- Användning av minst fyra olika spädningsfaktorer för de enstaka replikaten av ett prov

magellan kontrollerar hela layouten och väljer alla funna spädningsserier.

# 11.5.2 Curve Parameter Calculation (Beräkning av kurvparameter)

Kurvparametrar beräknas för varje funnen spädningsserie med hjälp av algoritmen **Four Parameters – Marquardt (Fyra parametrar – Marquardt)**. Om felet uppstår görs beräkningen om med hjälp av algoritmen **Four Parameters**. Se kapitel 11.4 Typer av analys av standardkurvor för mer detaljerad information om de algoritmer som nämns ovan.



# 11.5.3 Calculation of IC Values (Beräkning av IC-värden)

Först väljs den maximala intensiteten  $(I_{max})$  och den lägsta intensiteten  $(I_{min})$  i varje spädningsserie.

# Calculation of IC-intensity (Beräkning av IC-intensitet)

Om 0 %-värdet är inställt på intensiteten "Intensitet 0"

$$I_{IC} = \frac{ICx}{100.0} * I_{\max}$$

Om 0 %-värdet är inställt på "Min. intensitet för spädningsserie"

$$I_{IC} = I_{\min} + \frac{(I_{\max} - I_{\min}) * ICx}{100.0}$$

# Beräkning av spädning vid ICx

Med hjälp av de kurvparametrar som beräknats i steg 12.6.2 och den beräknade IC-intensiteten beräknas den resulterande spädningen.

*dilution* = 
$$C * (\frac{A - I_{IC}}{I_{IC} - D})^{\frac{1}{B}}$$

# 12. Exempel på tillämpning

# 12.1 Steg-för-steg exempel: Kvantitativ ELISA

Ett steg-för-steg-exempel (kvantitativt test) på hur man skapar en metod i **magellan** finns i det här kapitlet. Genom att följa instruktionerna lär du dig att definiera utvärderingar från en provsatsbeskrivning i **magellan**.

Metoden skapades med en Sunrise i demoläge, men det är också möjligt att definiera metoderna med ett annat Tecan-instrument anslutet, om det kan mäta absorbans.

# 12.1.1 Provsatsbeskrivning

I tillverkarens provsatsbeskrivning av kvantitativa IgM – Antikroppsdetektioner – ELISA finns följande instruktioner: Plattlayout

	1	2	3	4	5	6	7	8	9	10	11	12
A	BLK	C3	S1									
В	NC	C4	S2									
С	NC	C4	S2									
D	C1	C5	S3									
Е	C1	C5	S3									
F	C2	C6										
G	C2	C6										
Н	C3	S1										

BLK = Blank, NC = Negativ kontroll, C1–C6 = Kalibratorer (Standarder),

S1–S... = Prov

# Mätning och utvärdering

Läs av plattan vid en våglängd på 492 nm, referens vid 620 nm.

Blankreducera plattan med brunn A1.

Kalibratorers koncentration (Standarder):

Kalibrator 1	5 UA/ml
Kalibrator 2	10 UA/ml
Kalibrator 3	20 UA/ml
Kalibrator 4	40 UA/ml
Kalibrator 5	80 UA/ml
Kalibrator 6	160 UA/ml

Efter reduktionen för blanka ritas de optiska densiteterna (OD 492–OD620) upp mot respektive koncentration. Den regressionslinje som går genom dessa punkter utgör standardkurvan.



Tolkning av testresultaten:

IgM < 18 UA/ml	Negativ
18 UA/ml <= IgM < 22 UA/ml	Gränsfall
IgM >= 22 UA/ml	Positiv

Den beräknade IgM-koncentrationen hos de båda negativa kontrollerna måste understiga 8 UA/ml.

# Datahantering

Efter mätningen lagras datafilen (arbetsytan) automatiskt och en rapport med mätparametrar, plattlayout, blanka värden, standardkurva, IgM-koncentrationer, definition av cutoff, kvalitativa resultat för proverna och valideringar skapas.

Dessutom lagras layouten och de kvalitativa resultaten som ASCII-filer.

# 12.1.2 Create a Method (Skapa en metod)

I dialogrutan Wizard List (Guide-listan), välj Create/edit a method (Skapa/redigera en metod) och klicka på OK. Klicka på Next (Nästa) på Välkomstsidan på guiden Create/edit a method (Skapa/redigera en metod) och dialogrutan Select a File (Välj en fil) visas. Välj Create new (Skapa ny).

# **Measurement Parameters (Mätparameterar)**

Klicka på Nästa (Nästa) och dialogrutan Measurement parameters (Mätparametrar) visas.

Jreate/Edit a Method	>
Measur	rement type
	Endpoint measurement
	O Kinetic measurement
	Well kinetic measurement
	O Multilabel measurement
	Measurement parameters
- /	
Help	
Cancel <<< Back	

Välj **Endpoint measurement (Slutpunktsmätning)** och klicka på knappen **Measurement parameters (Mätparametrar)**. Följande dialogruta visas:



easurem	nent Paramete	r (SUNRISE Sim	ulation)		×
General	Wavelengths	Meas. Params	Temperature	Shaking	
۲	) Fixed wavelen	gth			
C	) Absorbance so	can	Set full wavelen	gth range	7.5
Wave	elength	2			
Me Ref	asurement wav	elength: ngth:	492 620	nm	
				OK Ca	incel Help

På fliken General (Allmänt) väljer du Absorbance (Absorbans).

På fliken Wavelengths (Våglängder) väljer du 492 nm som mätningsvåglängd och 620 nm som referensvåglängd.

Klicka på OK för att återgå till dialogrutan Measurement parameter (Mätparameter).

Klicka på **Define evaluation (Definiera utvärdering)** och fönstret **Plate layout (Plattlayout)** visas.





# Definiera layout

Definiera plattlayouten med hjälp av dialogrutan Well Assignment (Tilldelning av brunn) på höger sida av skärmen.

I grupprutan Identifiers (Identifierare) välj BL (Blank).

I rutan Experimental group (Försöksgrupp) kvarstår numret 1.

Välj All (Alla) i grupprutan Replicates (Replikat).



Klicka på väl A1 som sedan markeras med en röd kant.

Klicka på **Fill selection (Fyll urval)** och brunnen märks med den utvalda identifieringstypen.



Anmärkning En enstaka brunn kan också fyllas genom att dubbelklicka på den.

Välj nu följande inställningar i dialogrutan **Well Assignment (Tilldelning av brunn)**:

I grupprutan Identifiers (Identifierare) välj NC (Negativ kontroll)

I rutan Experimental group (Försöksgrupp) kvarstår numret 1.

I grupprutan Replicates (Replikat), väljs All (Alla).

Börja vid brunn **B1**, klicka och dra musen till **C1**. Brunnarna **B1** till **C1** markeras sedan med en röd kant.

Klicka på **Fill selection (Fyll urval)** och brunnarna märks med den utvalda identifieringstypen.

Därefter måste kalibratorer (standarder) tilldelas brunnarna **D1** till **G2**. Välj följande inställningar i dialogrutan **Well Assignment (Tilldelning av brunn)**:

I grupprutan Identifiers (Identifierare) välj ST (Standard).

I rutan Experimental group (Försöksgrupp) kvarstår numret 1.

l grupprutan Replicates (Replikat), välj mellan Fix number (Fast nummer) och All (Alla).

Fix number (Fast nummer)



Aktiveras endast för standarder och prover där ID:n kan användas. Om knappen **Fix number (Fast nummer)** är aktiv kan ett nummer skrivas in i motsvarande textfält. Detta nummer anger hur många replikat som är avsedda för denna metod. I de valda brunnarna skapas det angivna antalet replikat för varje ID. Därför måste antalet valda brunnar vara en multipel av det angivna antalet replikat.

#### All (Alla)

Alla valda brunnar definieras som replikat. Om ett befintligt ID-nummer för proverna och standarderna väljs, läggs de utvalda brunnarna till som replikat till de befintliga replikaten. Med alla andra identifieringstyper läggs de utvalda brunnarna till som replikat till de befintliga replikaten.

Två pilknappar anger riktningen för replikat och ID-nummersekvensens riktning (horisontell eller vertikal).

I det här exemplet, välj Fix Number (Fast nummer) och 2.

I rutan ID-Number (ID-nummer) och i grupprutan Replicates (Replikat) välj vertical arrows (vertikala pilar).

Välj sedan brunnarna D1 till G2 och klicka på Fill selection (Fyll urval).



Anmärkning Välj brunnarna enligt följande: Börja vid brunn D1, klicka och dra musen över de önskade brunnarna till H1. Håll sedan ner kontrolltangenten (Ctrl) och dra musen över de önskade brunnarna från A2 till G2.

Plate Layout (Plattlayouten) ser ut på följande sätt:



Klicka **Select all unused (Välj alla oanvända)** från verktygsfältet för att välja alla tomma brunnar på plattan. Håll sedan ner kontrolltangenten (Ctrl) och klicka på brunnen **H12** så att den förblir blank och omarkerad.

I dialogrutan Well Assignment (Tilldelning av brunn), välj SM (Sample) (Prov) under Identifiers (Identifierare).



I rutan Experimental group (Försöksgrupp) kvarstår numret 1.

I grupprutan Replicates (Replikat), välj Fix number (Fast nummer) och 2.

l rutan ID-Number (ID-nummer), lämnar du det som 1 och i grupprutan Replicates (Replikat), välj vertical arrows (vertikala pilar). Klicka sedan på Fill selection (Fyll urval). Layoutdefinitionen är slutförd.

# Transformationer

I kontrollfältet till vänster i fönstret väljer du nästa alternativ, Add new transformation... (Lägg till ny transformation) från elementet Transformed data (Transformerad data), för att definiera blankreducering.



En dialogruta kommer att fråga dig om du vill definiera en blankreducering. Klicka på **Yes (Ja)**.

Välj **Difference Data (Differensdata)** i rutan **Input data (Indata).** Som standard heter transformationen Transformation1 (se kontrollfältet). Om du tidigare har bekräftat definitionen av en blankreducering namnger programmet den automatiskt **Blank reduction (Blankreducering)**.

l rutan **Formula (Formel)** visas **x-BL1** automatiskt för denna blankreducering, där x refererar till det aktuella indatavärdet i en brunn och där BL1 är medelvärdet av de blanka brunnarna i försöksgrupp 1.

Välj nu hela plattan genom att klicka på -tecknet i mikroplattans övre vänstra hörn och klicka på den gröna bocken bredvid formelfönstret för att tilldela transformationerna till brunnarna. För ytterligare detaljer och förklaringar om definitionen och tilldelningen av transformationer, se kapitel 4.3.5 Transformed data (Transformerad data): Lägg till ny transformation. Följande fönster visas:

kent data (Utranere	e data		(v)									~ [A	alatile daria (hulfigile	data artic) 🔹
frederi		1	2	3	4	5	6	7	8	9	10	11	12	
- (2) 	A	8L1 1/1 x-8L1	5T1_3 22 x-BL1	SM1_1 2/2 x-BL1	SM1_5 2/2 x-BL1	SM1_9 2/2 x-BL1	SM1_13 2/2 x-BL1	SM1_17 2/2 x-BL1	SM1_21 2/2 x-BL1	SM1_25 2/2 x-BL1	SM1_29 2/2 x-BL1	5M1_33 2/2 x-8L1	SM1_37 2/2 x-BL1	
		NC1 1/2 x-BL1	ST1_4 1/2 x-BL1	SM1_2 1/2 x-BL1	SM1_6 1/2 x-BL1	SM1_10 1/2 x-BL1	SM1_14 1/2 x-BL1	SM1_18 1/2 x-BL1	SM1_22 1/2 x-BL1	SM1_26 1/2 x-BL1	SM1_30 1/2 x-BL1	SM1_34 1/2 x-8L1	SM1_38 1/2 x-BL1	
i handing		NC1 2/2 x-BL1	ST1_4 2/2 x-BL1	SM1_2 2/2 x-BL1	SM1_6 2/2 x-BL1	SM1_10 2/2 x-BL1	SM1_14 2/2 x-BL1	SM1_18 2/2 x-BL1	SM1_22 2/2 x-BL1	SM1_26 2/2 x-BL1	SM1_30 2/2 x-8L1	SM1_34 2/2 x-8L1	SM1_38 2/2 x-BL1	
		571_1 1/2 x-BL1	ST1_5 1/2 x-BL1	SM1_3 1/2 x-BL1	SM1_7 1/2 x-BL1	SM1_11 1/2 x-BL1	SM1_15 1/2 x-BL1	SM1_19 1/2 x-BL1	SM1_23 1/2 x-BL1	SM1_27 1/2 x-BL1	SM1_31 1/2 x-BL1	SM1_35 1/2 x-8L1	5M1_39 1/2 x-BL1	
		ST1_1 2/2 x-BL1	ST1_5 2/2 x-BL1	SM1_3 2/2 x-BL1	SM1_7 2/2 x-BL1	SM1_11 2/2 x-BL1	SM1_15 2/2 x-BL1	5M1_19 2/2 x-BL1	SM1_23 2/2 x-BL1	SM1_27 2/2 x-BL1	SM1_31 2/2 x-BL1	SM1_35 2/2 x-BL1	SM1_39 2/2 x-BL1	
	F	ST1_2 1/2 x-BL1	ST1_6 1/2 x-BL1	SM1_4 1/2 x-BL1	SM1_8 1/2 x-BL1	SM1_12 1/2 x-BL1	SM1_16 1/2 x-BL1	SM1_20 1/2 x-BL1	SM1_24 1/2 x-BL1	SM1_28 1/2 x-BL1	SM1_32 1/2 x-BL1	SM1_36 1/2 x-8L1	SM1_40 1/2 x-BL1	
		ST1_2 2/2 x-BL1	ST1_6 272 x-BL1	5M1_4 2/2 x-BL1	SM1_8 2/2 x-BL1	SM1_12 2/2 x-BL1	SM1_10 272 x-BL1	SM1_20 2/2 x-BL1	SM1_24 2/2 x-BL1	SM1_28 2/2 x-BL1	SM1_32 2/2 x-BL1	SM1_30 272 x-BL1	SM1_40 2/2 x-BL1	AN Austria
		ST1_3 1/2 x-BL1	SM1_1 1/2 x-BL1	SM1_5 1/2 x-BL1	SM1_9 1/2 x-8L1	SM1_13 1/2 x-BL1	SM1_17 1/2 x-BL1	SM1_21 1/2 x-BL1	SM1_25 1/2 x-BL1	SM1_29 1/2 x-BL1	SM1_33 1/2 x-BL1	SM1_37 1/2 x-8L1		TEC



SM1_9	Prov, försöksgrupp nummer 1, prov-ID nummer 4.
2/2	Antalet replikat är 2, det totala antalet replikat är 2.
x-BL1 eller 1.000	Tilldelad transformation x-BL1 (när Transformation är vald) eller spädningsfaktor 1 (när Conc., Dil,. Refvalues är valda). (Koncentration/spädning/referensvärde)

I varje brunn visas följande information (exempel brunn A5):

# Concentration / Dilution / Reference Value Definition (Definition av koncentration/spädning/referensvärde)

I kontrollfältet väljer du **Conc., Dil., Ref.-value (Konc., Dil., Ref.-värde)** från elementet **Metodlayout** för att definiera respektive värden enligt beskrivningen i testutrustningen.

Kalibrator 1	5 UA/ml
Kalibrator 2	10 UA/ml
Kalibrator 3	20 UA/ml
Kalibrator 4	40 UA/ml
Kalibrator 5	80 UA/ml
Kalibrator 6	160 UA/ml

Se till att ST väljs i listan Select Identifier (Välj identifierare).

I listan **Identifier (Identifierare)**, visas en lista av standarderna Grupp 1 visas. I den motsvarande **Concentration (Koncentration)**-rutan **ST1\_1** skriver du in siffran 5, och i rutan **Unit (Enhet)** skriver du in UA/ml. I den motsvarande **Concentration (Koncentration)**-rutan **ST1\_2** skriver du in siffran 10. Enheten behöver bara definieras en gång och gäller då för alla standarder. Skriv in värdena för ST1\_3 till ST1\_6 på samma sätt.

Skärmen som visar plattlayouten och koncentrationen visas:

Create/Edit a Method														×
Undo Parte NVL 1001. Zoon														
Select bardier     Select b	0	1	2	3	4	5	6	7	8	9	10	11	12	
Bok reductor         Image: Construction in the image: Construction in	A	BL1 1/1 1	ST1_3 2/2 20	SM1_1 2/2 1	SM1_5 2/2 1	SM1_9 2/2 1	SM1_13 2/2 1	SM1_17 2/2 1	SM1_21 2/2 1	SM1_25 2/2 1	SM1_29 2/2 1	SM1_33 2/2 1	SM1_37 2/2 1	
Conc. transforment d	B	NC1 1/2 1		SM1_2 1/2 1	SM1_6 1/2 1	SM1_10 1/2 1	SM1_14 1/2 1	SM1_18 1/2 1	SM1_22 1/2 1	SM1_26 1/2 1	SM1_30 1/2 1	SM1_34 1/2 1	SM1_38 1/2 1	
CC Nellinan Core landing Data bandhag Data coot Annual da bandhag Annual da bandhag	C	NC1 2/2 1	ST1_4 2/2 40	SM1_2 2/2 1	SM1_6 2/2 1	SM1_10 2/2 1	SM1_14 2/2 1	SM1_18 2/2 1	SM1_22 2/2 1	SM1_26 2/2 1	SM1_30 2/2 1	SM1_34 2/2 1	SM1_38 2/2 1	
Versional and a second and a se	D	ST1_1 1/2 5	ST1_5 1/2 80	SM1_3 1/2 1	SM1_7 1/2 1	SM1_11 1/2 1	SM1_15 1/2 1	SM1_19 1/2 1	SM1_23 1/2 1	SM1_27 1/2 1	SM1_31 1/2 1	SM1_35 1/2 1	SM1_39 1/2 1	
Urat [UAVAL]		9T1_1 2/2 5	ST1_5 2/2 80	SM1_3 2/2 1	SM1_7 2/2 1	SM1_11 2/2 1	SM1_15 2/2 1	SM1_19 2/2 1	SM1_23 2/2 1	SM1_27 2/2 1	SM1_31 2/2 1	SM1_35 2/2 1	SM1_39 2/2 1	
	F	ST1_2 1/2 10	ST1_6 1/2 160	SM1_4 1/2 1	SM1_8 1/2 1	SM1_12 1/2 1	SM1_16 1/2 1	SM1_20 1/2 1	SM1_24 1/2 1	SM1_28 1/2 1	SM1_32 1/2 1	SM1_36 1/2 1	SM1_40 1/2 1	.e
Aust	G	ST1_2 2/2 10	ST1_6 2/2 160	SM1_4 2/2 1	SM1_8 2/2 1	SM1_12 2/2 1	SM1_16 2/2 1	SM1_20 2/2 1	SM1_24 2/2 1	SM1_28 2/2 1	SM1_32 2/2 1	SM1_36 2/2 1	SM1_40 2/2 1	CAN Austr
	H	511_3 1/2 20	3M1_1 1/2 1	зм1_5 1/2 1	SM1_9 1/2 1	1/2 1/2	1/2 1/2	1/2 1/2	1/2 1/2	1/2 1/2	5M1_33 1/2 1	5MT_37 1/2 1		ΤĒ
The Carol Core Las													NEX	



# Standardkurva

l kontrollfältet, klicka på **Standard curve (Standardkurva)** från elementet **Concentrations (Koncentrationer)** för att definiera lämplig standardkurva.

Följande är i provsatsbeskrivning:

Efter reduktionen för blanka ritas de optiska densiteterna (OD 492–OD 620) upp mot respektive koncentration. Den regressionslinje som går genom dessa punkter utgör standardkurvan.

Välj Blank reduction (Blankreducering) som indata på fliken Data.

Standard curve	
Data Analysis type	Intercepts Axis Graph
Input data:	Blank reduction
Standards I	irom Layout
🔵 Standards I	rom ext. file: Select
🔘 Standards I	rom experimental group:
🔿 No Standar	d curve graph
Additional	concentrations

På fliken Analysis type (Analystyp), välj linjär regression.

	Analysis Type	
	O Point to point	
	Linear regression	Determine
	O Non-linear regression	
	O Cubic spline	endulend()
	O Akima	
	O Polynomial	
	O Four parameters	
	O Four parameters Marquardt	More
	O Five parameters	
	O LogitLog	
[	Include (0,0)	
Γ	Extrapolation factor: 1	



På fliken Axis (Axel), definierar du märkningen och skalningen av axlarna enligt nedan:

a Analysis type Intercepts Axis	Graph
X-axis	
Label: Concentration [UA/m]	11
Color:	Log. Scaling
Auto select range	
ORange	Min.: Max.:
🗹 Grid	Color: Style:
Y-axis	
Label: Blank reduction	
Color:	Log. Scaling
<ul> <li>Auto select range</li> </ul>	
⊖ Range	Min.: Max.:
🗹 Grid	Color: Style:

På fliken **Graph (Diagram)** kan du göra inställningar för diagramrubrik, kurvor, teckensnitt samt diagramvisning.

Analysis type Intercepts Axis	Graph		
Title			
Label: IgM ELISA			
Color:			
Curves			
	Label: Grp. 1		
Color:	Hide curve		
Symbol:	Line width:		
Size: Medium 🗸			
Font	Display		
Small	🖂 Legend	Intercepts	
O Medium	🗹 Base points	Error bars	



# Define Cutoffs (Definiera cutoffs)

Välj **Cutoff definition (Definition av cutoff)** i kontrollfältet från elementet**Evaluate data (Utvärdera data)** för att definiera gränserna för den kvalitativa utvärderingen.

Provsatsbeskrivningen innehåller följande instruktioner:

Tolkning av testresultaten:

IgM < 18 UA/mI	Negativ
18 UA/ml <= IgM < 22 UA/ml	Gränsfall
IgM >= 22 UA/mI	Positiv

Använd följande procedur för att definiera lämpliga cutoffs:

I rutan Ingångsdata väljer du Medelkoncentration. (UA/m).

Tabellen **Cutoffs** representerar en skala som visar höga respektive låga änden för **Limits (Gränser)** och **Labels (Benämningar)**. I **Limits (Gränser)**, skriv in 22 som det första (högre) gränsvärdet och 18 som det andra (lägre) gränsvärdet.

I Labels (Benämningar), skriver du in de olika testtolkningarna (Positive (Positiv), Intermediate (Gränsfall) och Negative (Negativ) i respektive ruta. Använd färgpaletten för att tilldela en färg:

Positiv – Röd

Gränsfall- Blå

Negativ - Grön

Skärmen innehåller följande:

Input data:	Mean conc. (UA/r	nL) 🗸	
Cutoffs			
Colors	Labels	Limits	
	Positive		
	Intermedia	22	
	▼ Negative	18	
		·	
	↓	·	
Formula inpu	ut		
Variable	C	Iperators Functions	
RI 1	~	t v and v	Cutoff results selection

Klicka på **Cutoff results selection (Urval Cutoff-resultat)** för att välja de identifieringstyper för vilka Cutoff-resultaten ska visas.


### Define QC Validations (Definiera QC-valideringar)

Klicka på **QC Validations (QC-valideringar)** i kontrollfältet från elementet **Evaluate data (Utvärdera data)**. Kriterierna för valideringen definieras så att giltigheten hos testresultaten kan garanteras.

I detta exempel måste följande krav vara uppfyllda:

Den beräknade IgM-koncentrationen hos båda negativa kontroller måste vara mindre än 8 UA/ml.

I rutan Input data (Indata) välj Single conc. (Enstaka koncentration) (UA/ml).

På den första raden skriver du **NC1\_1<8** eller ange formeln med hjälp av de tillgängliga **variabler**, **operatorer** och **funktioner**.



Anmärkning NC1\_1 betyder negativ kontroll för försöksgrupp 1, replikat 1.

På andra raden skriver du in NC1\_2<8.

Dialogrutan QC Validations (QC-valideringar) visas enligt följande:

3324	Input data:	Single conc. (U/	A/mL)	~	Validation group: 1	
Т	Validation	Conditions				1
1	NC1_1<8					
2	NC1_2<8					
3						
4						
5						
6	-					
7						
8						



### 12. Exempel på tillämpning

#### Organisera utskriven rapport

Klicka på **Printed report (Utskriven rapport)** i kontrollfältet från elementet **Data handling (Datahantering)**. Följande skärm visas:

Print as O List	You can use drag & drop to create a report. Pick an item from the available data list and drag it into the selected data list. Drop it into the appropriate position inside the report.
Available data:	Selected data:
Instrument data   Reduced data   Difference data - Mean   Difference data - Standard de   Difference data - Variation co   Transformed data   Transformed data   Oncentrations   Uaulitative Results   Image: Nethod layout   Method layout   UC Validation criteria   Measurement parameters   R Remarks   Fror protocol	Insert Append Ap
E_ Audit Trail E_ Signature	

På fliken **Data selection (Dataurval)**, finns alla data som är tillgängliga för rapporten i rutan **Available data (Tillgängliga data)**. Med hjälp av knapparna **Insert and Append (Infoga och bifoga)**, kan data överföras till rutan **Selected data (Utvalda data)**. Du kan även flytta data genom att klicka och dra med musen.

I rutan **Print as (Skriv ut som)** kan du välja mellan att skriva ut data i form av en matris eller som en lista med en specifik orientering.

I det här exemplet bör en rapport med mätparametrar, plattlayout, blanka värden, standardkurva, IgM-koncentrationer, definition av cutoff, kvalitativa resultat av proverna och valideringar skapas.

Innan rapporten skapas, ska standardinställningen Matrix Difference data (Differensdata i matris) tas bort från rutan Selected data (Utvalda data) rutan. Därefter återstår endast Measurement parameters (Mätparameterar) i rutan Selected data (Utvalda data).

Välj Method layout (Metodlayout)/Layout i rutan Available data (Tillgängliga data) och bifoga den som en matris till rapporten genom att klicka på Append (Bifoga). Infoga därefter Blank reduction (Blankreducering), Mean conc. (Medelkoncentration) (UA/ml) samt Cutoff results (Cutoff-resultat) i denna matris genom att välja de motsvarande elementen och sedan klicka på Insert (Infoga).



Append Graph (Bifoga diagram): Standardkurva, definition av cutoff och Kriterier för QC-validering till de utvalda data. Därmed är dataurvalet och förberedelsedelen av rapportdefinitionen slutförd. Dialogrutan Printed Report (Utskriven rapport) ska se ut på följande sätt:



På flikarna **Header (Sidhuvud)** och **Footer (Sidfot)** definierar du layouten för rapportens sidhuvud och sidfot (se kapitel 4.3.13 Datahantering: Utskriven rapport för mer information).



#### Data Export (Dataexport)

Klicka på **Data export (Dataexport)** i kontrollfältet från elementet **Data handling** (**Datahantering**). I detta exempel ska layout samt de cutoff-resultaten sparas som en ASCII-fil. Välj Layout och Cutoff-resultaten från fönstret Available data (Tillgängliga data). Klicka på  $\rightarrow$  pilen för att infoga dem i fönstret Selected data (Utvalda data). Skärmen visar följande information:

⊕ 🛍 Instrument data 🔨 🔨		Layout	Up
		Lutoff results	Down
			Dow
Difference data - Variation coefficier			
Transformed data			
	->		
Cutoff results	<-		
Sample IDs			
Method layout			
Strip method names			
Original Concentrations			
Dilution factors			
			tan Tan
11	200000		



Anmärkning Data som exporteras bör alltid innehålla layout eller prov-ID.

### Automated Data Handling (Automatiserad datahantering)

l kontrollfältet väljer du Automated Data Handling (Automatiserad datahantering) från elementet Data handling (Datahantering).

Automatic	
load sample ID list	More
z save workspace	More
export to ASCII file	
export to ASTM file (LIS)	More
export to Sample Tracking	
export to Excel	
_ print	
-	

Välj export to ASCII file (exportera till ASCII-fil), och view results after measurements (visa resultat efter mätningar). I magellan Tracker är Save Workspace (Spara arbetsyta) vald som standard och kan inte ändras.





## Spara metoden

Klicka på **Next (Nästa)** för att öppna fönstret **Save as (Spara som)**. Ange ett filnamn för metoden och fyll i övriga fält vid behov.

Create/Edit a Method	×						
Save in: mth Filename: Application Example_en.mt File remarks: Audit trail comment: Organize favorites Audit	Name       Remarks       Status         Application Example_ch       Unsigned         Application Example_en       Unsigned         Application Example_en       Unsigned         Application Example_fr       Unsigned         Application Example_fi       Unsigned         Application Example_fi       Unsigned         Application Example_fi       Unsigned         Application Example_po       Unsigned         Application Example_po       Unsigned         Application Example_ru       Unsigned         Application Example_po       Unsigned         Mathitabel_SUNRISE       Unsigned         Multilabel_SUNRISE       Unsigned         Multilabel_SUNRISE       Unsigned         Spectrum_SUNRISE       Unsigned         Ital       Signatures						
Cancel Cancel Cancel	Run this method now SAVE&FINISH						
Filename (Filnamn) Fältet File remarks (Filkommentarer)	De kommentarer som anges här kommer att sparas och visas tillsammans med filnamnet.						
Textfältet Audit trail comment (Audit trail kommentar)	De kommentarer som anges här kommer att sparas i Audit trail. Detta alternativ är bara tillgängligt i <b>magellan</b> <b>Tracker</b> .						
Knappen Audit trail	Dialogrutan <b>Audit Trail</b> visas. Detta alternativ är bara tillgängligt i <b>magellan Tracker</b> .						
Knappen Organize Favorites (Organisera favoriter)	ialogrutan <b>Organize Favorites (Organisera favoriter)</b> sas. Se kapitel 6.4 Start Favorite (Starta Favorit) – Organize avorites (Organisera favoriter)).						
Knappen Signatures (Underskrifter)	Dialogrutan <b>Signature (Underskrift)</b> visas. Detta alternativ är bara tillgängligt i <b>magellan Tracker</b> .						
Method password (Lösenord för metod)	Ange ett lösenord för metoden om du vill skydda den metod som ska sparas (se kapitel 4.4.1 Lösenordsskydd av metoder).						
Kryssrutan Run this method now (Kör denna	Metoden körs omedelbart efter att du har avslutat guiden.						

metoden nu)



#### Audit Trail

Klicka på Audit Trail för att visa dialogrutan Audit Trail:

Administrator (Ad         2018-03-06         09:02:44         Created           Administrator (Ad         2018-03-06         09:13:59         Modified	

Audit trail visar en lista över alla ändringar som gjorts i metoden. Varje inmatning består av användaren (namn och fullständigt namn), datum och tid för ändringen, om filen skapades eller ändrades och eventuella kommentarer till audit trail.

Klicka på **Print Preview... (Förhandsgranskning av utskrift)** för att få en förhandsvisning av filen. För att jämföra en metod med tidigare versioner måste en utskrift göras, eftersom två fönster för förhandsgranskning av utskrift inte kan öppnas samtidigt.

## 12.1.3 Kör metoden

Om du har valt **Run this method now (Kör denna metod nu)** i dialogrutan **Save** as (Spara som) i Create/edit a method wizard (Guiden Skapa/redigera en metod), kommer dialogrutan **Start Measurement Wizard/Start Measurement** (Starta mätning) att visas när du klickar på **Save (Spara)**.

t Measurement					
Measurement			Measurement parameters		
Workspace:	06032018-001.wsp	Arb. cycle kin	Measurement mode: Absorbance		
Method: Application Example_en		Modify layout Measurement wavelength: 492 nm Beference wavelength: 620 nm			
Sample ID list:		Insert	Read mode: Normal		
Instrument	acker	Current: 32.5 °C			
Plate in	Temp. control	Target: 37.0 °C			
Movements	Injector control	Optimize Z-position			
Please note:		-/			
	/				
Help					
Cancel	<<< Back		START V		



Klicka på **Start** för att starta mätningen. Nu skapas automatiskt en arbetsyta som innehåller all tidigare angiven information, och som kommer att samla upp alla mätvärden. Under tiden som mätningen pågår visas en statusruta som indikerar hur mätningen fortskrider.

När mätningen har slutförts visas dialogrutan **Results (Resultat)**, i vilken alla resultat och beräkningar visas.

## 12.1.4 Evaluate the Result (Utvärdera resultat)

**Evaluate results (Utvärdera resultaten)** gör det möjligt för dig att visa och utvärdera rådata. Du kan se utvärderingsparametrarna och utvärdera data på nytt.

Det här avsnittet guidar dig genom guiden **Evaluate results (Utvärdera resultat)** med hjälp av den exempelarbetsytan som skapats genom att köra den kvantitativa ELISA-metoden.



#### Anmärkning

Exempelfiler visas automatiskt i metodlistan i magellan Standard. Dessa filer finns tillgängliga för magellan Tracker i standarddatasökvägen för data och måste konverteras.

Klicka på Evaluate results (Utvärdera resultat) genom dialogrutan Wizard List (Guide-lista).

När du klickar på **Next** (Nästa) på sidan **Welcome** (Välkomstsidan) av guiden **Evaluate Results wizard (Guiden Utvärdera resultat)** visas dialogrutan **Select a file (Välj en fil)**.

Välj arbetsytan **Quantitative Elisa example\_Sunrise.wsp** i fillistan och klicka på **Make your selection (Gör ditt val)**. Beräkningarna utförs och följande fönstret plattlayout visas:

50. <b>26</b>		1	2	3	4	5	6	7	8	9	10	11		
· · · · · · · · · · · · · · · · · · ·	A	BL1 1/1	ST1_3 2/2	SM1_1 2/2	SM1_5 2/2	SM1_9 2/2	SM1_13 2/2	SM1_17 2/2	SM1_21 2/2	SM1_25 2/2	SM1_29 2/2	SM1_33 2/2	SM1_37 2/2	
the native		NC1 1/2	ST1_4 1/2	SM1_2 1/2	SM1_6 1/2	SM1_10 1/2	SM1_14 1/2	SM1_18 1/2	SM1_22 1/2	SM1_26 1/2	SM1_30 1/2	SM1_34 1/2	SM1_38 1/2	
2	C	NC1 2/2	ST1_4 2/2	SM1_2 2/2	SM1_6 2/2	SM1_10 2/2	SM1_14 2/2	SM1_18 2/2	SM1_22 2/2	SM1_26 2/2	SM1_30 2/2	SM1_34 2/2	SM1_38 2/2	
2		ST1_1 1/2	0.646 ST1_5 1/2	0.128 SM1_3 1/2 0.287	0.289 SM1_7 1/2	-0.001 SM1_11 1/2	0.123 SM1_15 1/2	-0.001 SM1_19 1/2	0.312 SM1_23 1/2 0.497	0.129 SM1_27 1/2	-0.002 SM1_31 1/2	-0.001 SM1_35 1/2	0.327 SM1_30 1/2	
28		ST1_1 2/2	ST1_5 2/2	SM1_3 2/2 0.282	SM1_7 2/2 0.501	SM1_11 2/2 0.499	5M1_15 2/2	SM1_19 2/2	SM1_23 2/2	SM1_27 2/2	SM1_31 2/2 0.201	SM1_35 2/2 0.385	SM1_39 2/2 0.301	
		ST1_2 1/2	ST1_6 1/2	SM1_4 1/2	SM1_8 1/2	SM1_12 1/2	SM1_10 1/2	SM1_20 1/2	SM1_24 1/2	SM1_28 1/2	SM1_32 1/2	SM1_36 1/2	SM1_40 1/2	
54		ST1_2 2/2	ST1_6 2/2 2.345	SM1_4 2/2	SM1_8 2/2 0.270	SM1_12 2/2 0.303	SM1_16 2/2	SM1_20 2/2 0.775	SM1_24 2/2	SM1_28 2/2	SM1_32 2/2	SM1_36 2/2 0.402	SM1_40 2/2 0.41	4 Austria
	H	ST1_3 1/2	SM1_1 1/2	SM1_5 1/2	SM1_9 1/2	SM1_13 1/2 0.512	SM1_17 1/2	SM1_21 1/2	SM1_25 1/2	SM1_29 1/2	SM1_33 1/2	SM1_37 1/2		TECA
													-	
Measurement parameters  SUMRES Instrument secial number 1212004177 Measurement model: Absolvance Measurement unovidingtis: 482 nm Reference ausedingtis: 482 nm														

I varje enstaka brunn visas det beräknade värdet. Fönstret plattlayouten ändras på motsvarande sätt, beroende på det valda elementet i kontrollfältet. Parametrar och inställningar kan ändras med hjälp av elementen i kontrollfältet. Om metoden ska ändras, klicka på fliken **Edit method (Redigera metod)**.



Klicka i brunnen med höger musknapp och en sammanhangsberoende meny visas:

SM1_1	SM1_5	SM1_9	SM				
2/2	2/2	2/2	2				
0.293	Summary						
SM1_	Details						
1/2	Edit						
0.119	Edit kinetic settings						
CM1	Copy kinetic settings						
	Paste kinetic settings						
2/2	Graph: Multilabel						
0.128	Graph: Kinetics						
SM1	Graph: Multilabel kinetics						
4/0	Graph: FLT curves						
1/2	Graph: Spectra						
0.287	Graph: Dilution	n series					
SM1_	Mask/Unmask	selection					
2/2	Show/Hide lay	vout					

Om du väljer **Summary (Sammanfattning)** visas följande fönster med detaljerad information om definitionen och inställningarna för den valda brunnen:

- 🔁 🚺 (	Concentrations	1
	Single conc. (UA/mL)	
	17.121	
	Average single conc. (UA/mL)	
	🖥 Mean conc. (UA/mL)	
. 🗄 🚺 🤅	Itatistics	
]	👠 s - Measurement data	
	🛕 v - Measurement data	
	👠 s - Reference data	
	🛕 v - Reference data	
	👠 s - Difference data	
	🛕 v - Difference data	
	🛕 s - Blank reduction	
	🛕 v - Blank reduction	
	👠 s - Average single conc. (UA/mL)	
	🛕 v - Average single conc. (UA/mL)	
	Up	
	Down	

Klicka på **Finish (Slutför)** i fönstret plattlayout och **Save as (Spara som)** visas, där du kan ange ett filnamn och kommentarer. Klicka på den lilla **Save (Spara)**knappen till vänster i fönstret för att spara filen. Du kan fortsätta arbeta med metoden eller arbetsytan. Klicka på knappen **Save (Spara)** till höger längst ner på skärmen för att spara filen och stänga guiden. Programmet går tillbaka till guide-listan.



## 12.1.5 Sammanfattning av definitionen av kvantitativ ELISA i magellan

#### 1. Dra bort blanka värden

#### Definitioner i magellan

Klicka på Add new transformation (Lägg till ny transformation) i kontrollfältet och ett fönster kommer att fråga dig om du vill definiera en Blank reduction (Blankreducering). Klicka på Yes (Ja) och formeln Blank reduction (Blankreducering) tilldelas automatiskt alla brunnar.

#### 2. Definiera koncentrationer

Definitioner i magellan (Kontrollfält – Method layout/ Conc.-, Dil.-, Ref.values) (Metodlayout/Koncentration/spädning/referensvärde)

Utvald identifierare: ST Enhet: UA/ml

5	(ST1 1Standard 1 första försöksgrupp)
10	(ST1_2Standard 2 första försöksgrupp)
20	(ST1_3Standard 3 första försöksgrupp)
40	(ST1_4Standard 4 första försöksgrupp)
80	(ST1_5Standard 5 första försöksgrupp)
160	(ST1_6Standard 6 första försöksgrupp)
	5 10 20 40 80 160

#### 3. Definiera standardkurva

Definitioner i magellan (Kontrollfält – Concentrations/ Standard curve) (Koncentrationer/Standardkurva)

Indata	blankreducering
Analystyp	linjär regression
X-axel	linjär
Y-axel	linjär

### 4. Define Cutoffs (Definiera cutoffs)

Definitioner i magellan (Kontrollfält – Evaluate data/ Cutoff definition) (Utvärdera data/Definition av cutoff)

• Indata: Mean conc. (UA/ml)

Gränser 22

18

Positiv>=22>intermediär>=18>negativ Icke-kompetitivt test#

## 5. QC Validation (QC-validering)

Definitioner i magellan (Kontrollfält – Evaluate data/ Cutoff definition) (Utvärdera data/Definition av cutoff)

• Indata: Single conc. (UA/ml)

Valideringsvillkor 1 NC1\_1<8 Valideringsvillkor 2 NC1\_2<8

NC1\_1.....Negativ kontroll första replikat första försöksgruppen NC1\_2.....Negativ kontroll andra replikat första försöksgruppen

# 13. Ordlista med termer

Term	Definition
Average single conc. (???)	Koncentrationen beräknad genom att använda medelvärdet av de enstaka koncentrationerna.
Basis OD	Kinetisk parameter: Bas OD-värde för beräkningen av starttiden
Basis OD %	Kinetisk parameter: Bas OD-värde för beräkningen av starttiden i %
Urklipp	Urklipp är det medium genom vilket program i Windows utbyter data med varandra. Data kan väljas i ett Windows-program genom att klippa eller kopiera den till urklipp och sedan läggas till i ett annat program genom att klistra in den från urklipp.
Korrelationskoefficient	Korrelationskoefficienten anger styrkan och riktningen av ett linjärt samband mellan två slumpmässiga variabler.
Cutoff-gräns	Med hjälp av Cutoff-gräns kan användaren definiera gränserna mellan två tillstånd (t.ex. positivt eller gränsfall). Dessa kriterier används vid utvärderingen av resultaten.
Cutoff-resultat	Namnet på det cutoff-intervallet som värdet ligger inom enligt definitionen under definition av definition
Definiton av cutoff	Definition av alla cutoff-intervall och beräknade gränser
Spädningsfaktorer	Definierade spädningsfaktorer för proverna och kontrollerna. En spädningsfaktor på 2 motsvarar en spädning 1:2.
Spädningsserie	Ett prov med replikat i olika spädningar.
Anpassningsgrad	1 minus de genomsnittliga relativa kvadratiska avvikelserna för baspunkterna från kurvan.
Diagram	Diagram kan visas för kinetiska mätningar, enzymkinetik, mätningar med multipla inmärkningar, spädningsserier eller standardkurvor.
Dold	Visas om brunnsdata döljs vid utskrift
HUID	Hardware Unit Identification Number (Identifikationsnummer för maskinvara)
IC50	Den spädning/koncentration som resulterar i 50 % av det maximala responsen
Ogiltig	Värdet är ogiltigt, ingen beräkning möjlig
Graph: Kinetic	Diagram över kinetiska mätningar
Lamp Low	Inga värden från mätningen på grund av fel i absorbansinstrumentet.
Layout, Plattlayout	Definierar var prover eller kontroller placeras på mikroplattan.
Max. slope OD/hr	Kinetisk parameter: Maximal lutning av de kinetiska kurvorna per timme.
Max. slope OD /min	Kinetisk parameter: Maximal lutning av de kinetiska kurvorna per minut.
Max. slope OD /sec	Kinetisk parameter: Maximal lutning av de kinetiska kurvorna per sekund.
Maximum OD	Kinetisk parameter: Maximala värdet för de kinetiska kurvorna
Mean slope OD /hr	Kinetisk parameter: Genomsnittlig lutning av kinetiska kurvorna per timme

#### 13. Ordlista med termer



Term	Definition
Mean slope OD /min	Kinetisk parameter: Genomsnittlig lutning av kinetiska kurvorna per minut.
Mean slope OD /sec	Kinetisk parameter: Genomsnittlig lutning av kinetiska kurvorna per sekund.
Mean. conc (???)	Koncentration beräknad från medelvärdet av replikat av indata.
Mätdata	Absorbansmätning med dubbla våglängder: Data som mäts med hjälp av mätfiltret
Mätparameterar	Definierar mätläge, våglängd, plattstorlek, skakning osv.
Mättyp	Mättypen kan vara slutpunktsmätning, kinetisk mätning, mätningar med multipla inmärkningar eller kinetisk mätning i brunn.
Method (Metod)	Metoderna innehåller mätparametrarna och utvärderingsdefinitionen. När du kör en metod öppnas en arbetsyta som innehåller de uppmätta och beräknade data.
Minimum OD	Kinetisk parameter: Minsta värde för de kinetiska kurvorna
MultPt	Standardkurvan är inte monoton och ger mer än en koncentration vid givna indata.
NoCalc	Inget värde har returnerats från beräkningen.
Ursprungliga koncentrationer	Koncentrationer i den standardkurva som definieras i metoden.
Pipetteringsstatus	Om en prov-ID-lista importeras från ett pipetteringsprogram kan pipetteringsstatusen visas.
QC-validering	Kriterierna för QC-validering definieras i metoden och anger om en mätning är giltig eller ogiltig. Kriterierna kan till exempel ange om mätvärdena ligger för långt ifrån varandra eller om de avviker för mycket från ett förväntat värde. Programmet varnar automatiskt användaren om kriterierna inte uppfylls.
Rådata	Data som mäts av instrumentet.
Referensdata	Absorbansmätning med dubbla våglängder: Data som uppmätts med referensfiltret
Resultatstatistik	Sammanfattning av antalet värden i de olika cutoff-intervallerna.
Sample ID List (Prov-ID-lista)	Prov-ID:n tilldelas varje brunn på grund av att den tillhörande sonden kan identifieras. ID:n är vanligtvis streckkoder som importeras från prov-ID-listor som lagras av ett pipetteringprogram.
Prov-ID:n	ID:n för proverna
Single. conc (???)	Koncentration beräknad från indata för varje replikat
Test	I tidigare versioner innehöll ett test utvärderingsinställningarna men inte mätparametrarna. Tester stöds inte längre i <b>magellan</b> och ersätts av mer kraftfulla metoder.
Time Basis OD	Kinetisk parameter: Tid till dess att basvärdet har uppnåtts
Time Basis OD	Kinetisk parameter: Tid tills basvärdet % har uppnåtts
Time Basis to Onset OD	Kinetisk parameter: Tid mellan basvärde och startvärde
Time Basis to Onset OD %	Kinetisk parameter: Tid mellan basvärdet och startvärdet %
Time max. slope sec	Kinetisk parameter: Kinetisk parameter: Tidpunkt för den maximala lutningen
Time maximum OD	Kinetisk parameter: Tid till dess att det maximala värdet har uppnåtts

Term	Definition
Time minimum OD	Kinetisk parameter: Tid till dess att miniminivån är uppnådd
Time Onset OD	Kinetisk parameter: Tid till dess att det definierade startvärdet har uppnåtts
Time Onset OD %	Kinetisk parameter: Tid till dess att den definierade starttiden i % har uppnåtts.
Tidpunkter	Tidsstämplar för de enstaka mätningarna i en kinetisk mätning.
Ej tillgänglig	De begärda data är inte tillgängliga.
Oanvänd	Inga data uppmätt; ingen layout definierad i denna brunn.
Användarprompt	Användarprompter är referenser som ställs in i metoden. De visas före mätningen och användaren måste lägga till text i dem. Denna text kommer sedan att ingå i en utskrift.
Valideringskriterier	Sammanfattning av resultaten av valideringsvillkoren
Brunnsposition	Brunnens namn, till exempel: A1, A2, …
Workspace (Arbetsyta)	Alla tillgängliga data i programmet finns i arbetsytan, t.ex. mätdata, utskriftsdefinitionen och metoddefinitionen. Arbetsytan används för att ladda in metoder och utföra mätningar.
!	Föregår värden som har eliminerats under beräkningen
#	Föregår koncentrationsvärden som har beräknats med hjälp av extrapolering och som ligger utanför standardkurvans intervall
()	Parenteser omger värden som har dolts.
*	En asterisk markerar värden som har uppmätts med alternativet "Use gain regulation" (Använd förstärkningsreglering), som korrigerar (= sänker) förstärkningen.
??? – Mean	Beräknade medelvärden (till exempel: Rådata – medelvärde)
<pre>??? - Standard deviation s - ???</pre>	Beräknade standardavvikelser (till exempel Rådata – Standardavvikelse eller s – Rådata)
<pre>??? - Variation coefficient v - ???</pre>	Beräknade variationskoefficienter (till exempel: Rådata – Variationskoefficient eller v – Rådata)
~	Läggs till värden som har redigerats eller simulerats.
<min< th=""><th>Den beräknade koncentrationen är lägre än det minsta</th></min<>	Den beräknade koncentrationen är lägre än det minsta
>Max	Den beräknade koncentrationen överskrider det maximala
<blank></blank>	Utskriven rapport: Infoga en tom matris eller tabellutskrift
<page break=""></page>	Utskriven rapport: Skriv ut nästa element på nästa sida
<separator></separator>	Utskriven rapport: Skriv ut en linje mellan två element
X	Symbol <b>x</b> hänvisar till det aktuella värdet i en brunn.
concX	Symbolen <b>concX</b> avser koncentrationen av standarden i den aktuella brunnen.
dilX	Symbolen <b>dilX</b> avser spädningen av provet eller kontrollen i den aktuella brunnen.
'???'!	Tillgänglig dataset om fler än ett set indata ska användas för beräkningarna.



#### 13. Ordlista med termer

Term	Definition
[]	Index anger de olika cyklerna i en kinetisk mätning, medan [0] anger den första cykeln.
*	Asterisken markerar identifierare som har sätts som alias.



# Register

## A

A	
abs(argument)	187
Akima	199
Alternativ for E-postmeddelande	168
Alternativ for kopiering/export	156
Ä	
Ändra lösenord	25
Α	
Ansluta ett instrument	
Använd fördefinierad metod	109, 111
Användaradministration	
Alternativ	166
Användaradministration (magellan Tracker)	
Ändra användare	170
Inloggning	170
Användargränssnitt	17
Användarprompt	
Användarrättigheter	172
Användarrättigheter (magellan Standard)	171
Arbiträr kinetikcykel	113
arean under den kinetiska kurvan	61
Arkivfiler	148
ASTM filexport	
ASTM-fil	83, 85
Audit Trail	177
Audit trail för användaradministration	165
Autofyll urval	
Automatiserad datahantering	
Automatisk arkivering	149
avg(argument)	188

## В

Beräkna på nytt med en annan metod	128
Beräkningar	
Reducering av spektradata	182
Beräkningar	181
Beräkningar	181
Bifoga underskrift	145

# С

calcAlways(argument)	193
countDeleted(arg1; arg2)	
Cutoff	
Cybersäkerhet	

## D

Dataexport7	71
Datahantering7	71
Definiera en ny identifierare4	18
Definiera en plattlayout4	45
Definiera filterhållare	34
Definiera konstanter5	59
Definiera utvärdering4	12
Definition av ASTM-avgränsare	33
Diagram för enzymkinetik 13	31
Diagram för standardkurva 13	35
Diagram med multipla inmärkningar 13	38
Dialogrutan Spektra	41
Dialogrutan Välkommen2	23
Driftskvalificering OQ	16

## Е

eliminate (arg1; arg2; arg3; arg4) eliminatePerc (arg1; arg2; arg3; arg4) eliminateRange (arg1; arg2; arg3; arg4) eller – logiskt uttryck	190 190 191 186
Enzymkinetik	63
Erhålla rådata	. 109, 110
Exempel på kvantitativ ELISA	
Exempel på tillämpningar	
exp(argument)	
Export av ASCII-fil	
Export av ASTM-fil	, 123, 124
Export av ASTM-fil (LIS)	
Export av ASTM-filer	
Export av Excel	
Export av metod	
Exportera	
till ASCII	74
till Excel	75
Exportera till LIS	

## F

Fem parametrar – Marquardt	203
Filhantening	10 / ۱
Fliken Kopiera	156
Fliken Övrigt	150 150
Fliken Plattwy	159 158
Fliken Sökvägar	150
Förster plattlavout	120
Fönstret plattlavout	123 43
Förflyttning	
nlatthållare filterhållare	33
Förhandsgranskning av utskrift	
Formel	183
Andra funktioner	193
Fliminationsfunktioner	190
Funktioner	186
Grundläggande funktioner	187
Hur man skriver en formel	183
Logiska uttryck	
Spektrafunktioner	193
Statistiska funktioner	188
Variabler	183
frac(argument)	187
Fyra parametrar	202
Fyra parametrar – Marquardt	203
• •	

# G

Genvägslista Godkännande	
Granskning	
Guide	
Standardelement	22
Guide-lista	17
Guiden Registrera	
Guiden Skapa/redigera en prov-ID-lista	91
н	
HUID	14, 28, 178
I	
ICx-beräkning identifierare	52 48, 49

	400
if() then() else() – logiska uttryck	
ignore()	193
Ikonen Ovrigt	147
Alternativ	154
Ikoner	18
Importera en prov-ID-lista	99
Importera rådata	150
Importera/redigera en prov-ID-lista	93
In(argument)	187
Infoga en prov-ID-lista	115
Information om magellan	176
inloganing	170
Inloggning	25
Inloggningar	-
misslvckade	
Installation	
Programvara	
Installationskyalificering	13
Inställningar för ASTM-export	
Inställningar för plattvy	158
Inställningar för prov-ID-lista	Q1
Instrumentkentrell	
isinvalid()	193

## Κ

Kinetiska diagram	139
Kinetiska mätningar	41
Kinetiska parametrar	131
Kinetiska transformationer	63, 68
Klistra in från ASCII-format	128
Klistra in i ASCII-format	
Knappen Hjälp	22
Kompatibilitet med läsare	12
Komponenter och termer	19
Koncentration, spädning och referensvärden	50
Koncentrationer	132
Konvertera dokument	151
Konvertera från	152
Konvertera till	151
Kopiera i ASCII-format	128
Kopiera till Excel	128
korrelationskoefficient	.203, 204
Kriterier för filval	120
Kubisk spline	199
Kvalitativa resultat	136

## L

Lägg till HUID	178
Lägg till ny transformation	56
Lägg till/ändra användare (magellan Standard) .	170
Lägg till/ändra användare (magellan Tracker)	162
Lägg till/ändra roll	164
Lås program	166
Levenberg-Marquardt-algoritmen	203
Levenberg-Marquardt-metoden	203
Levy-Jennings-diagram	125
lg(argument)	187
Licensiera magellan	28
Log(argument)	187
Logiska uttryck	
eller	186
if() then() else()	186
och	186
Lösenord	
Alternativ	167
Lutning	59
-	



# Μ

magellan Standard	12
magellan Tracker	12
Mapphantering	20
Mätning med multipla inmärkningar för absorbans	40
Mätningar med multipla inmärkningar	38
Mätparameterar	37
Mätparametrar	37
Mätstatus1	16
Mättyper - SUNRISE	37
max(argument)1	89
maxAvg(argument)1	89
mean(argument)1	88
median(argument)1	88
medianPlate()1	88
Menyer i fliken Utvärdera resultat 1	23
Metodanteckningar	86
Metoder med flera plattor	88
Metodlayout1	36
min(argument)1	89
Min./Max	61
minAvg(argument)1	89
Misslyckade inloggningar 1	66

## Ν

Nummerformat 86
-----------------

#### 0

186
227
111
121
136

## Ρ

Plattlavout	45
PointwiseCV(argument)	
Programmet låst	25
Programvara	
Avinstallation	
Installation	
Prov-ID	136
Prov-ID-lista	
importera	

# Q

QC platta-till-platta	71,	125
QC-validering	70,	136

## R

Redigera användarprompt	112
Redigera brunn	
Redigera konstanter	112
Reducerade data	130
Reducering av kinetiska data	59
Reducering av spektradata	
Registreringsformulär	30
round(argument)	187

## S

Sammanfattning av användaradministrationen 168 Sammanfattning av brunn
Detaljer 137
Dialogrutan Diagram-kinetik 139
Dialogrutan Diagram-multipla inmärkningar 138
Dialogrutan Diagram-Spektra 141



Dialogrutan Redigera	138
Sammanfattning av en brunn	43, 137
Särskilda karaktärer	130
Signera en fil	145
Skapa/redigera en metod	35
Skriv ut	124
Skrivarinställningar	124
Slutpunktsmätningar	40
SMTP-server	168
Spädningsserie	135
Spara de utvärderade resultaten	143
Spara metoden	
Spara prov-ID-listan	107
Spektrum för en skanningsmätning	141
sqr(argument)	188
sqrt(argument)	188
Standardelement	
Standardkurva	63, 132
Standardsökvägar	155
Starta Favorit	109
Starta magellan	24
Starta mätning	109, 112
Starttider	61
stddev(argument)	189
Sum(argument)	189
Symboler för filtyper	20
System Audit Trail	179
Systemåterställning	14

## Т

Teckensnitt för utskrift Temperaturkontroll Testa e-post Tilldela alias Tillgängliga data Tillverkare	
Transformerad data	131
Typer av analys av standardkurvor	196

### U

Underskrift	145
Godkännande	
Granskning	
Ursprungliga lösenord	25
Utskriven rapport	76
Utvärdera data	68
Utvärdera resultat	119, 143

#### **V** Viktoir

Viktning	
Anpassning med fyra/fem parametrar –	
Marquardt/polynomanpassning	204
W	
Westgard®-regler	27

# Varumärken

Följande produktnamn och eventuella registrerade eller oregistrerade varumärken som nämns i denna handbok används endast som referens och förblir respektive ägares exklusiva egendom:

- [Produktnamn]<sup>TM</sup> (infinite®)
   Tecan® och Tecan-logotypen är registrerade varumärken som tillhör Tecan Group Ltd., Männedorf, Schweiz
- DNA Expert<sup>TM</sup> är ett registrerat varumärke som tillhör Techcomp Ltd., Hong Kong, Kina
- Agilent® är ett registrerat varumärke som tillhör Agilent Technologies, Inc., Santa Clara, CA, USA
- AIR LIQUIDE<sup>™</sup> är ett registrerat varumärke som tillhör AIR LIQUIDE, S.A., Paris, Frankrike
- Aseptisol® är ett registrerat varumärke som tillhör BODE Chemie GmbH & Co. KG, Hamburg, Tyskland
- Bacillol® är ett registrerat varumärke som tillhör BODE Chemie GmbH & Co. KG, Hamburg, Tyskland
- Costar®, Corning® och NBS<sup>TM</sup> är registrerade varumärken som tillhör Corning Incorporated, Corning, NY, USA
- Greiner®, µClear®, Lumitrac<sup>™</sup> och Fluotrac<sup>™</sup> är registrerade varumärken som tillhör Greiner Labortechnik GmbH, Frickenhausen, Tyskland
- HTRF® är ett registrerat varumärke som tillhör Cisbio International, Frankrike
- Invitrogen<sup>TM</sup> är ett registrerat varumärke som tillhör Invitrogen Corporation, Carlsbad, CA, USA
- Lysetol® och Gigasept® (tidigare Lysetol) är registrerade varumärken som tillhör Schülke & Mayr GmbH, Norderstedt, Tyskland
- Microcide® är ett registrerat varumärke som tillhör Global Biotechnologies Inc., Portland, Maine, USA
- Microman® är ett registrerat varumärke som tillhör Gilson, Inc., Middleton, WI, USA
- Pentium® är ett registrerat varumärke som tillhör Intel Corporation, Santa Clara, CA, USA
- Invitrogen® och PanVera® är registrerade varumärken som tillhör Invitrogen Corporation Carlsbad, CA, USA
- Windows®, MS DOS®, Visual Basic® och Excel® är registrerade varumärken som tillhör Microsoft Corporation, Redmond, WA, USA
- Hamamatsu® är ett registrerat varumärke som tillhör HAMAMATSU Photonics K.K. [IR], Hamamatsu City, Japan
- NUNC<sup>TM</sup> och Matrix är registrerade varumärken som tillhör Thermo Fisher Scientific, Waltham, MA, USA
- Polyfiltronics® är ett registrerat varumärke som tillhör Whatman International Ltd.
- Dynex® är ett registrerat varumärke som tillhör magellan Biosciences, Chelmsford, MA, USA
- Labsystem® är ett registrerat varumärke som tillhör Labsystem kft., Budapest, Ungern
- BRET<sup>2</sup>®, DeepBlueC® och PerkinElmer® är registrerade varumärken som tillhör PerkinElmer, Inc., Waltham, Massachusetts, USA
- Chroma-Glo<sup>TM</sup> är ett registrerat varumärke som tillhör Promega Corporation, Madison, WI 53711 USA
- MycoAlert® är ett registrerat varumärke som tillhör Cambrex Corporation, East Rutherford, NJ, USA

# **Tecan kundsupport**

Kontakta Tecans lokala kundsupportorganisation om du har några frågor eller behöver teknisk support för din Tecan-produkt. Kontaktuppgifter finns på <u>http://www.tecan.com/</u>.

Innan du kontaktar Tecan för att få produktsupport ber vi dig ha följande information tillgänglig för att vi ska kunna hjälpa dig på bästa möjliga sätt (se typskylten):

- Produktens modellnamn
- Produktens serienummer
- Programvara och programversion (om tillämpligt)
- Beskrivning av problemet och kontaktperson
- Datum och klockslag då problemet inträffade
- Steg som redan vidtagits för att korrigera problemet
- Dina kontaktuppgifter (telefonnummer, e-postadress, faxnummer osv.).



# Licensavtal för slutanvändare av programvara

VIKTIGT - LÄS NOGGRANT: Detta Licensavtal ("LICENSAVTAL") är ett bindande avtal mellan Er (fysisk eller juridisk person) och Tecan Österrrike GmbH ("Tecan") för den "proprietary" programvaruprodukten som identifieras ovan, vilken inkluderar programvara ("PROGRAMVARA") och kan inkludera tillhörande media, skriftligt material, "online"- eller elektronisk dokumentation ("DOKUMENTATION"). Denna PROGRAMVARA ställs till Ert förfogande endast under de villkor som framgår av detta LICENSAVTAL. Genom att installera, kopiera eller på annat sätt använda PROGRAMVARAN (vilket som först inträffar), har Ni accepterat att följa villkoren i detta LICENSAVTAL. Om Ni inte accepterar villkoren i detta LICENSAVTAL, har Ni inte rätt att använda PROGRAMVARAN.

#### PROGRAMVARULICENS

PROGRAMVARAN och DOKUMENTATIONEN är skyddade av upphovsrättslig lagstiftning och andra immaterialrättsliga lagar och konventioner. PROGRAMVARAN säljs inte, den licensieras.

#### 1.1.LICENSMEDGIVANDE

Detta LICENSAVTAL medger en icke-exklusiv och icke-överförbar rätt att använda PROGRAMVARAN i enlighet med instruktionerna och procedurerna i DOKUMENTATIONEN enbart för Ert interna bruk. Ni får installera och använda det antal kopior av PROGRAMVARAN, på de antal datorer och arbetsstationer anslutna till det antal enheter, som specificerats i DOKUMENTATIONEN till detta LICENSAVTAL i enlighet med proceduren för programvaruinstallation som beskrivs i DOKUMENTATIONEN.

Ni har rätt att antingen göra en kopia av PROGRAMVARAN enbart för arkivändamål till stöd för Er användning av PROGRAMVARAN på en enstaka dator, eller överföra PROGRAMVARAN till en enstaka hårddisk förutsatt att Ni behåller originalet enbart för säkerhets- eller arkivändamål. Ni har inte rätt att reproducera eller distribuera PROGRAMVARAN på något annat sätt, inklusive utlåning eller leasing.

PROGRAMVARAN skall endast användas tillsammans med de enheter som specificeras i dokumentationen och får inte användas med några andra enheter.

#### 2. LICENSAVGIFT

Licensavgiften för rätten att använda PROGRAMVARAN fastställs i DOKUMENTATIONEN.

#### 3. BESKRIVNING AV ANDRA RÄTTIGHETER OCH BEGRÄNSNINGAR

Ni har inte rätt att utföra Reverse Engineering, dekompilering eller disassemblering av PROGRAMVARAN, med undantag av och endast i den utsträckning att en sådan aktivitet uttryckligen är tillåten enligt lag.

PROGRAMVARAN licensieras som en enstaka produkt. Dess komponenter får inte delas upp för användning på fler än en dator eller arbetsstation.

Tecan kan tillhandahålla supporttjänster relaterade till PROGRAMVARAN ("SUPPORTTJÄNSTER"). Användning av SUPPORTTJÄNSTER styrs av Tecan's policy och program vilka beskrivs i användarhandboken, i "online"-DOKUMENTATIONEN och/eller i annat material som tillhandahålls av Tecan. Alla tilläggsprogram och all programkod som överlämnas till Er som en del av SUPPORTTJÄNSTERNA ska betraktas som en del av PROGRAMVARAN och underkastas villkoren i detta LICENSAVTAL. Beträffande den den tekniska information som Ni överlämnar till Tecan som en del av SUPPORTTJÄNSTER, kan Tecan använda sådan information för sina affärssyften, inklusive produktsupport och utveckling. Tecan kommer inte att utnyttja denna tekniska information på ett sådant sätt att det kan identifiera Er personligen.

Ni har inte rätt att överföra någon av Era rättigheter som faller under detta LICENSAVTAL utan ett föregående skriftligt medgivande från Tecan.

Oaktat varje annan rättighet, kan Tecan upphäva detta LICENSAVTAL om Ni underlåter att efterfölja villkoren i LICENSAVTALET. I sådant fall har Ni inte längre rätt att använda PROGRAMVARAN. Ni är då skyldig att förstöra alla kopior av PROGRAMVARAN, DOKUMENTATIONEN och alla dess komponenter inom fjorton dagar.

#### 4. UPPHOVSRÄTT

Alla äganderätter och upphovsrätter i och till PROGRAMVARAN (inklusive, men inte begränsat till, de diagram, bilder, fotografier, animationer, videor, ljud, musik, text och "applets" som införlivats i PROGRAMVARAN), DOKUMENTATIONEN och alla kopior av PROGRAMVARAN och DOKUMENTATIONEN ägs av Tecan eller dess underleverantörer och är licensierade till Tecan. PROGRAMVARAN och DOKUMENTATIONEN är skyddade av upphovsrättslig lagstiftning och andra immaterialrättsliga lagar. Därför måste Ni behandla PROGRAMVARAN och DOKUMENTATIONEN på samma sätt som varje annat upphovsrättsskyddat material.

#### 5. VILLKOR FÖR ANVÄNDNING

Användningen av PROGRAMVARAN skall ske under åtlydande av följande villkor:

- PROGRAMVARAN skall enbart användas tillsammans med enheter som specificerats i DOKUMENTATIONEN;
- enbart den senaste versionen av PROGRAMVARAN skall användas;
- PROGRAMVARAN får inte modifieras eller förorsakas att bli modifierad.

Det åligger Er personligen att själv beakta alla myndighetskrav angående användandet av denna PROGRAMVARA.



#### 6. FÖRSIKTIGHETSMÅTT VID ANVÄNDNING

Innan PROGRAMVARAN blev färdig produkt, undergick PROGRAMVARAN en serie av tester för att mäta tillförlitligheten i de uppnådda resultaten.

Emellertid, med hänsyn taget till det stora antalet möjliga tillämpningar för vilka PROGRAMVARAN kan användas, har det inte varit möjligt att utföra sagda tester i en verklig arbetssituation.

De avtalsslutande parterna är överens om att det i praktiken inte är möjligt att ta fram databehandlingsprogram som är 100 % lämpliga för alla tillämpningar.

Därför rekommenderar vi starkt att Ni verifierar och validerar erhållna resultat innan PROGRAMVARAN används i en verklig arbetssituation varje gång den ska användas i en ny tillämpning.

Skulle det uppstå några tvivel angående resultaten, eller i sådant fall att felaktiga resultat kan påvisas, uppmanas Ni att omedelbart kontakta Tecan.

För en perfekt användning av PROGRAMVARAN påminns Ni om att:

- Ni måste ha erforderliga kunskaper;

- Ni måste, på sedvanligt sätt, använda skrivskydd och i synnerhet säkerhetskopiering, för att skydda datafiler från oförutsedda händelser såsom radering, överskrivning, virusinfektioner osv.

#### 7. BEGRÄNSAD GARANTI

PROGRAMVARAN står under garanti under en period av 90 dagar efter leveransdatum.

Under denna period kommer Tecan att vidta skäliga ansträngningar för att leverera, så snabbt som möjligt och på det sätt som Tecan finner lämpligt, en uppdaterad version av programmet utan varje betydande defekt som påvisats.

De åtgärder som nämns ovan utgör omfattningen av den garanti som medges Er.

Påpekande om Java-stöd. PROGRAMVARAN kan innehålla stöd för program skrivna i Java Java-teknologin är inte feltolerant och har inte utformats, tillverkats eller avsetts för användning eller återförsäljning som en komponent i kontrollutrustning i farliga miljöer som kräver felsäker operation, som t ex styrsystem för kärnenergianläggningar, navigerings- eller kommunikationssystem för flygtrafik, flygledarsystem, direkt livsuppehållande maskiner eller vapensystem, i vilka en brist i Java-teknologin skulle kunna leda till döden, personskada eller svår fysisk eller miljömässig skada. Sun Microsystems, Inc. kräver att denna friskrivningsklausul inkluderas.

#### 8. BEGRÄNSAT ANSVAR

Tecan skall underkastas ett åtagande att utöva tillbörliga åtgärder.

Sagda åtagande gäller tillhandahållandet av programvaran i enlighet med den funktionalitet som beskrivits i dokumentationen av densamma, däremot ges ingen som helst garanti med hänsyn till de resultat som erhålls eller lämpligheten av PROGRAMVARAN för Ert ändamål.

Ni är ensam ansvarig för användningen av PROGRAMVARAN och för de resultat som erhålls. Ni har ett särskilt ansvar för att verifiera innehållet i resultaten och konsekvenserna därav, eftersom PROGRAMVARAN skall användas uteslutande under Ert överinseende, kontroll och ansvar.

TECAN SKALL INTE HÅLLAS ANSVARIG FÖR NÅGRA INDIREKTA ELLER PÅFÖLJANDE SKADOR ELLER EVENTUELL FÖRLORAD VINST, MINSKNING I INTÄKTER ELLER FÖRLUST AV DATA SOM ÅSAMKATS AV ER, ÄVEN OM TECAN BLIVIT INFORMERAT OM ATT NÄMNDA FÖRLUST ELLER SKADA KUNDE INTRÄFFA.

PROGRAMVARAN har undergått ett anti-virus-test. Dock skall inte Tecan hållas ansvarig om ett virus, som var omöjligt att upptäcka med hjälp av det anti-virus-verktyg som användes i den ovan nämnda testen, skulle visa sig vara närvarande.

SKULLE AV NÅGOT SOM HELST SKÄL TECAN HÅLLAS ANSVARIG INFÖR DOMSTOL, SKALL BELOPPET AV EVENTUELLT SKADESTÅND SOM TECAN ÅLÄGGS ATT UTBETALA, MED UNDANTAG FÖR UPPSÅTLIGT FEL ELLER GROV VÅRDSLÖSHET, INTE ÖVERSTIGA DEN SUMMA SOM MOTTAGITS AV TECAN I KRAFT AV DETTA LICENSAVTAL.

#### 9. KONFIDENTIALITET

Ni bekräftar härmed och accepterar att Ni inte kommer att delge PROGRAMVARAN eller DOKUMENTATIONEN till någon av Era anställda (med undantag av de som är ansvariga för programvaran) eller till någon tredje part utan föregående skriftligt medgivande från Tecan och att denna skyldighet till konfidentialitet överlever villkoren i detta LICENSAVTAL. Ni skall skydda PROGRAMVARAN och DOKUMENTATIONEN med största möjliga aktsamhet.

#### **10. UPPSÄGNING**

Om Ni underlåter att åtlyda någon av förpliktelserna i detta avtal, har Tecan rätten att när som helst säga upp LICENSAVTALET och omedelbart ta PROGRAMVARAN och DOKUMENTATIONEN och alla dess kopior, oavsett var de finns, i beslag utan föregående krav eller meddelande.

TECAN AUSTRIA GMBH, Untersbergstrasse 1a, A-5082 Grödig / Salzburg, Austria T +43 62 46 89 330, F +43 62 46 72 770, office.austria@tecan.com, www.tecan.com



# **Declaration of Conformity**

We, TECAN Austria GmbH herewith declare under our sole responsibility that the product identified as:

Product Type: Software

Model Designation: SW MAGELLAN V7.5

Article Numbers: 30143623, 30143624, 30143625, 30143626, 30143627, 30143628

Address: Tecan Austria GmbH Untersbergstr. 1A

A-5082 Grödig, Austria SRN: AT-MF-000020241

is in conformity with the provisions of the following EC Directive(s)/Regulation(s) when installed in accordance with the installation instructions contained in the product documentation:

#### **Regulation IVD-R**

on in vitro diagnostic devices

For products placed on the Swiss market the CH authorized representative is:

Tecan Schweiz AG Seestrasse 103, 8708 Männedorf, Switzerland



is in conformity with the relevant U.K. legislation for UKCA-marking when installed in accordance with the installation instructions contained in the product documentation:

#### Medical Devices Regulations 2002

Classification: Other device (all devices except Annex II and self-testing devices) Conformity assessment procedure: Annex III

For products placed on the United Kingdom market the **UK responsible person** is: Tecan UK Limited Theale Court, 11-13 High Street, Theale Reading, Berkshire, RG7 5AH United Kingdom

The current applicable versions of the directives and regulations as well as the list of applied standards which were taken in consideration can be found in separate CE & UK declarations of conformity.

These Instructions for Use and the included Declaration of Conformity are valid for all Magellan SW Packages with the article numbers listed above.