# **•**TECAN•

Användarhandbok för absorbansläsaren

# **INFINITE F50 PLUS**

och programvaran

# MAGELLAN



Dokumentnummer:	30225416
	2023-06
Dokumentutgåva:	1.2
Fast programvaruversion:	3.33
Programvaruversion Magellan:	7.5







#### VARNING

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

#### Observera

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 detta dokument.



#### 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 detta dokument 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 EU-godkännande

Se sista sidan i användarhandboken.

#### Om denna användarhandbok

Ursprungliga instruktioner. Detta dokument är avsedd som användarinstruktioner för absorbansläsaren INFINITE F50 PLUS, vilket är utformad för att mäta ljusabsorbansen (optisk densitet) på prover i mikroplattor med 96 brunnar. Det är avsett som referens och användarinstruktioner.

Detta dokument beskriver hur man:

- Installerar instrumentet
- Använder instrumentet
- Rengör och underhåller instrumentet

#### Anmärkning om skärmbilder

Det kan hända att versionsnumret som visas på skärmbilderna inte stämmer överens med numret på den aktuella utgåvan. Skärmdumpar 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, försiktighetsåtgärder och anmärkningar

Följande typer av meddelanden används i den här publikationen för att belysa viktig information eller för att varna användaren om eventuellt farliga situationer:



# Innehållsförteckning

1.	Säke	Säkerhet7		
	1.1	Instrum	entsäkerhet	7
2.	Allm	ant		
	2 1	hbaev&	användning	۰۹
	22	Funktio	nsprincin	
	2.3	Använd	arprofil	
		2.3.1	Professionell användare – administratörsnivå	
		2.3.2	Slutanvändare eller rutinanvändare	
		2.3.3	Servicetekniker	11
3.	Kom	nma igån	q	13
•	31	Unnack	aning och inspektion	13
	0.1	3.1.1	Inspektion av levererad förpackning	
		3.1.2	Steg för uppackning	
	3.2	Strömkr	rav	
	3.3	Miliökra		
	3.4	System	krav	15
	3.5	Slå PÅ i	instrumentet	16
	3.6	Progran	nvara	17
		3.6.1	Introduktion/avsedd användning	17
		3.6.2	Installationprocedur	17
		3.6.3	Installationskvalificering	18
		3.6.4	Börja arbeta med Magellan	19
	3.7	Magella	n - Measurement Parameter Editor (Redigerare för mätparametrar)	21
		3.7.1	Kontrollfält	22
		3.7.2	Workflow Pane (Panel för arbetsflöde)	26
		3.7.3	Infopanel	26
	3.8	Magella	n - Definiera mätningar	27
		3.8.1	Definiera slutpunktsmätningar	27
		3.8.2	Definiera mätningar med multipla inmärkningar	
		3.8.3	Definiera kinetiska mätningar	29
	~ ~	3.8.4	Dra in och slappa programelement	
	3.9	Optimer	ring for basta prestanda	32
		3.9.1	Tillyägagångagött	∠دعد مد
		3.9.2	Tillvagagangssatt	
4.	Inst	rumentfu	nktioner	33
	4.1	Instrum	entfunktioner	
		4.1.1	Skakning av mikroplatta	
	4.2	Instrum	entbeskrivning	
	4.3	Beskriv	ning av filterhjul	
	4.4	Instrum	entspecifikationer	<b>30</b>
		4.4.1	Aimanna specifikationer	0د 27
		4.4.2	Maispecifikationer	
	45	Instrum	enttillbehör	38
5	Kva	litotekont		20
5.	5 4	Interskom		
	5.1	Siälyko	ly ntroll	
	5.2	Driftelov	valificering (OQ)	
	5.5	531	Tester med MultiCheck	۵۵ مح
		5.3.2	Tester med mikroplattor	כם אריייייסא
		5.3.3	Instrumentprecision med flytande prover	
		5.3.4	Instrumentliniäritet med flytande prover	
6	Eve	mnel nå t		13
υ.			ווומווואָדוווואַ	43 40
	0.1 6.2	Stor för	19 r-stag avampal: Kvantitativ El ISA	43 43
	0.2	oleg-ioi	olog exemper. Itvanulativ LEIGA	



		6.2.1	Provsatsbeskrivning	43
		6.2.2	Create a Method (Skapa en metod)	45
		6.2.3	Organisera rapport	55
		6.2.4	Kör metoden	59
		6.2.5	Utvärdera resultat	60
		6.2.6	Sammanfattning av definitionen av kvantitativ ELISA i Magellan	62
7.	Ren	göring, ı	underhåll, och avfall	65
	7.1	Inledni	ng	65
	7.2	Rengö	ring av instrumentet	65
	7.3	Desinfe	ektion av instrument	66
		7.3.1	Desinfektionslösningar	66
		7.3.2	Desinfektionsprocedur	67
		7.3.3	Säkerhetsintyg	68
	7.4	Föreby	ggande underhållschema för INFINITE F50 PLUS	69
		7.4.1	Månatlig	69
		7.4.2	Vart fjärde år	69
	7.5	Filterby	yte och installation	69
		7.5.1	Steg för filterutbyte	70
		7.5.2	Definiera filter	73
	7.6	Avfall.		73
		7.6.1	Inledning	73
		7.6.2	Avfallshantering av förpackningsmaterial	73
		7.6.3	Avfallshantering av förbrukat material	73
		7.6.4	Avfallshantering av instrumentet	74
8.	Fels	ökning .		75
	8.1	Inledni	ng	75
		8.1.1	Tabell med felmeddelanden och felsökning	75
		8.1.2	Definition av 'Overflow (Överflöde)'	
		8.1.3	Strömavbrott	
9.	Förk	ortning	ar, varumärken och symboler	77
	9.1	Förkor	tningar	
	9.2	Varuma	ärken	
	9.3	Symbo	ler	
Reai	ster	-		81
Teca	n kun	dsuppo	rt	83



# 1. Säkerhet

# 1.1 Instrumentsäkerhet

- 1. Vidta alltid grundläggande försiktighetsåtgärder när du använder produkten för att minska risken för skada, brand eller elektriska stötar.
- 2. Läs och förstå all information i användarhandboken. Underlåtenhet att läsa, förstå och följa instruktionerna i detta dokument kan leda till att produkten skadas, att operatören skadas eller att instrumentet fungerar dåligt. Tecan ansvarar inte för skador som orsakats av felaktig hantering av enheten.
- 3. Observera alla VARNINGAR och uppmaningar till FÖRSIKTIGHET i detta dokument.
- 4. Koppla alltid bort enheten från huvudströmförsörjningen före rengöring och desinfektion.
- 5. Öppna aldrig instrumenthöljet.
- 6. Vidta lämpliga försiktighetsåtgärder för laboratorier, till exempel att bära skyddskläder (handskar, laboratorierock och skyddsglasögon) och tillämpning av godkända laboratorierutiner.



#### FÖRSIKTIGHET

OM INSTRUKTIONERNA I DENNA PUBLIKATION INTE FÖLJS KORREKT KAN INSTRUMENTET SKADAS. PROCEDURER KAN KANSKE INTE UTFÖRAS KORREKT OCH INSTRUMENTETS SÄKERHET KAN INTE GARANTERAS.

Det förutsätts att de som använder instrumentet, på grund av sin arbetslivserfarenhet, känner till nödvändiga försiktighetsåtgärder vid hantering av kemikalier och biologiska substanser.

Följ följande lagar och riktlinjer:

- Nationella arbetsskyddslagar
- Riktlinjer för att förhindra olyckor
- Säkerhetsdatablad från reagens-tillverkare

#### VARNING

BEROENDE PÅ TILLÄMPNINGARNA KAN DELAR AV INFINITE F50 PLUS HA KOMMIT I KONTAKT MED BIOLOGISKT/INFEKTERAT MATERIAL.

SE TILL ATT ENDAST KVALIFICERAD PERSONAL ANVÄNDER INSTRUMENTET. VID BEHOV AV SERVICE ELLER OM INSTRUMENTET FLYTTAS ELLER KASSERAS, SKA INSTRUMENTET ALLTID DESINFICERAS I ENLIGHET MED INSTRUKTIONERNA I DETTA DOKUMENT.

VIDTA LÄMPLIGA FÖRSIKTIGHETSÅTGÄRDER FÖR LABORATORIER, TILL EXEMPEL ATT BÄRA SKYDDSKLÄDER VID ARBETE MED POTENTIELLT SMITTFÖRANDE SUBSTANSER.







#### VARNING

INSTRUMENTET UPPFYLLER KRAVEN FÖR EMISSION OCH IMMUNITET SOM BESKRIVS I IEC 61326-2-6, MEN DEN ELEKTROMAGNETISKA MILJÖN BÖR UTVÄRDERAS INNAN INSTRUMENTET ANVÄNDS.

DET ÄR OPERATÖRENS ANSVAR ATT SE TILL ATT DEN ELEKTROMAGNETISKA MILJÖN ÄR LÄMPLIG FÖR INSTRUMENTET SÅ ATT INSTRUMENTET GER FÖRVÄNTADE RESULTAT.

ANVÄND INTE INSTRUMENTET I NÄRHETEN AV STARKA ELEKTROMAGNETISKA KÄLLOR (T.EX. OSKÄRMADE AVSIKTLIGA RF-KÄLLOR) EFTERSOM DETTA KAN STÖRA INSTRUMENTETS FUNKTION OCH LEDA TILL FELAKTIGA RESULTAT.



# 2. Allmänt

# 2.1 Avsedd användning

INFINITE F50 PLUS är en automatiserad absorbansläsare för mikroplattor med 96 brunnar, inklusive Magellan-programvara för professionell användning i ett laboratorium för mätning av ljusabsorption (optisk densitet) av homogena flytande medier för användning vid in vitro-diagnostik.

Instrumentet är avsett att användas främst vid in vitro-diagnostiska analyser av prover från människokroppen som levereras från en användarvald Enzyme-linked Assay (ELISA - Enzymkopplad immunadsorberande analys). Den specifika diagnostiska informationen och provtypen definieras av den valda analysen.

INFINITE F50 PLUS är avsedd för mätning och utvärdering av kvalitativa, semikvantitativa och kvantitativa analyser enligt planerade diagnostiska parametrar och instrumentspecifikationer.

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





Anmärkning Om INFINITE F50 PLUS absorbansläsare eller Magellan-programvaran ändras på något sätt är garantin inte längre giltig och instrumentet förlorar sitt regulatoriska efterföljande.

#### Anmärkning

Den verksamhetsansvariga får endast använda CE-märkta provsatser för tillämpningar inom klinisk diagnostik. Den verksamhetsansvariga måste säkerställa att kombinationen av en särskild CE-märkt provsats och den CE-märkta INFINITE F50 PLUS absorbansläsare och dess tillval har validerats att efterfölja IVD-direktivet 2017/746 eller andra relevanta nationella eller lokala regelverk.

Om INFINITE F50 PLUS absorbansläsare används på annat sätt än den "avsedda användning" som beskrivs ovan, eller om den används med annan programvara än Magellan, så uppfyller instrumentet inte längre IVD och användaren är ansvarig för användning och nödvändig validering.



Anmärkning

Resultat som erhållits från INFINITE F50 PLUS påverkas av korrekt användning av instrumentet och mikroplattor, i enlighet med instruktionerna i detta dokument, samt av de vätskor som används (reagenter, kemi). Användarhandboken, lagring och tillämpningar som inkluderar prover eller reagenter måste följas strikt. Resultaten måste därför tolkas noggrant.



Anmärkning Öppna aldrig instrumenthöljet, annars upphör garantin att gälla.



# 2.2 Funktionsprincip

INFINITE F50 PLUS absorbansläsare med åtta kanaler mäter den optiska densiteten (OD) av vätskor i mikroplattor med 96 brunnar.

Ljuset från LED-ljuskällan passerar genom ett optiskt filter i filterhjulet för att välja den våglängd som behövs för absorbansmätningen.

Efter passeringen genom filtret delas ljusstrålen upp i åtta optiska fibrer för att fokusera på åtta brunnar i mikroplattan. Ljusstrålarna som passerar genom mikroplattans brunnar fokuseras av åtta optiska linser och skickas sedan till åtta fotodioder för detektion.

Detta innebär att INFINITE F50 PLUS mäter åtta brunnar parallellt, vilket motsvarar en kolumn i en mikroplatta med 96 brunnar.

För att kontrollera och reglera LED-lampans ljusintensitet används ytterligare en optisk fiber (referenskanal), där ljuset passerar förbi mikroplattan.

De erhållna transmissionsvärdena omvandlas till värden för optisk densitet (OD) i enligt med följande formel:

Transmissi on T = 
$$\frac{I}{I_{\circ}}$$

I<sub>0</sub> = intensitet av incidentens ljus

 $I_0$  = intensitet av överförd ljus

OD är logaritmen av den omvända transmissionen.

$$OD = Log \frac{1}{T}$$



# 2.3 Användarprofil

## 2.3.1 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.

#### 2.3.2 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.

#### 2.3.3 Servicetekniker

Serviceteknikern är en person med lämplig teknisk utbildning och motsvarade färdigheter och erfarenhet. Om produkten behöver service eller underhåll kan personen identifiera och undvika farliga situationer.

Vana vid datorprogram och goda kunskaper i engelska krävs.



Anmärkning Information om utbildningsdatum, utbildningarnas längd och hur ofta de genomförs finns tillgänglig via kundsupport.

> Adress och telefonnummer finns på webben: http://www.tecan.com/customersupport



# 3. Komma igång

# 3.1 Uppackning och inspektion

#### 3.1.1 Inspektion av levererad förpackning

Det levererade instrumentet inkluderar:

- Extern strömförsörjning
- Elkabel
- USB-kabel f
   ör anslutning till extern dator
- Användarhandbok, PDF-filer på lagringsmedia
- USB-minne
  - Programvaran Magellan
  - o Användarhandbok
  - o Verktyg (t.ex. Adobe Reader)



Anmärkning För att undvika oönskade dataförluster eller angrepp av virus/skadlig attack får du aldrig ta bort skrivskyddet från USB-minnet.



#### FÖRSIKTIGHET

LÄSAREN HAR TESTATS MED DEN MEDFÖLJANDE USB-KABELN. OM EN ANNAN USB-KABEL ANVÄNDS KAN INSTRUMENTETS KORREKTA PRESTANDA INTE GARANTERAS.

## 3.1.2 Steg för uppackning

- 1. Inspektera förpackningen visuellt för skador innan den öppnas. Rapportera omedelbart eventuella skador.
- 2. Välj en plats att placera instrumentet på. Platsen bör vara plan, fri från vibrationer, skyddad mot direkt solljus samt fri från damm, lösningsmedel och sura ångor. Se till att det finns ett avstånd på minst 5 cm mellan instrumentet och väggen eller annan utrustning.
- 3. Lyft upp instrumentet ur kartongen och placera det på den valda platsen. Var försiktig när du lyfter upp instrumentet.
- 4. Inspektera instrumentet visuellt för eventuella lösa, böjda eller skadade delar. *Rapportera omedelbart eventuella skador.*
- 5. Jämför serienumret på instrumentets bakpanel med serienumret på följesedeln. Rapportera omedelbart eventuella avvikelser.
  - Rapponera officacioan eventacila avvikeiser.
- 6. Kontrollera instrumentets tillbehör och jämför med packsedeln.
- 7. Spara förpackningsmaterialet för eventuell senare frakt.



## 3.2 Strömkrav

Instrumentet känner automatiskt av den anslutna spänningen. Det är därför inte nödvändigt att göra några ändringar i spänningsområdet. Kontrollera spänningsspecifikationerna och se till att den spänning som levereras till instrumentet är korrekt enligt följande specifikationer:

#### Spänning:

Grundläggande instrumentet med AC-adapter:	100–240 V AC, 50/60 Hz
Grundläggande instrumentet utan AC-adapter:	24 V DC

Om den ovannämnda spänningen inte är tillgänglig i ditt land, kontakta Tecans lokala kundsupport.

Instrumentet får endast anslutas till strömkällor med skyddsjord.



#### FÖRSIKTIGHET ANVÄND INTE INSTRUMENTET I ETT FELAKTIGT

ANVAND INTE INSTRUMENTET I ETT FELAKTIGT SPÄNNINGSOMRÅDE. INSTRUMENTET SKADAS OM DEN SÄTTS PÅ MED FELAKTIG SPÄNNING.



#### FÖRSIKTIGHET

#### INSTRUMENTETS NÄTKABEL FÅR BARA ERSÄTTAS MED EN KABEL MED SAMMA SPECIFIKATIONER.

# 3.3 Miljökrav

Instrumentet bör placeras på en plan, slät yta som är fri från damm, lösningsmedel och sura ångor.

Vibrationer och direkt solljus skall undvikas för att säkerställa korrekta resultat.

#### Omgivningstemperatur:

Drift	15 °C till 35 °C (59 °F till 95 °F)
Förvaring	-30 °C till 60 °C (-22 °F till 140 °F)
Relativ luftfuktighet:	20-80 % icke-kondenserande vid driftstemperatur



# 3.4 Systemkrav

	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), 2019 (32-bitars), Microsoft Excel 365 (32-bitars) Endast 32-bitarsutgåvor stöds! Startutgåvor stöds INTE!	2010 (32-bitars) 2019 (32-bitars)



# 3.5 Slå PÅ instrumentet

De följande procedurer beskriver de nödvändiga stegen som krävs innan instrumentet sätts på.



#### FÖRSIKTIGHET

INNAN INSTRUMENTETS INSTALLATION OCH PÅSLAGNING SKA DET STÅ UTAN ATT VARA PÅSLAGET I TRE TIMMAR FÖR ATT ELIMINERA RISK FÖR KONDENSBILDNING SOM KAN ORSAKA KORTSLUTNING.

När de krav som nämns ovan har uppfyllts, utförs installationen enligt följande procedur:

- 1. Anslut instrumentet till den externa datorn med USB-kabeln.
- 2. Kontrollera att huvudströmbrytaren på instrumentets vänstra sida står i läge "OFF" (Av).
- 3. Sätt in elkabeln i eluttaget i den vänstra panelen.
- 4. Slå på instrumentet med huvudströmbrytaren på den vänstra panelen.

Instrumentet är klar att mäta mikroplattor efter installation av programvaran.



Anmärkning Kontrollera att mikroplattans position A1 är korrekt innan du påbörjar mätningar.

Mikroplattor kan endast mätas utan lock.

Stäng platthållarens skydd innan du påbörjar en mätning för att undvika att omgivande ljus påverkar resultaten.

Anmärkning Ta alltid bort mikroplattan från läsaren direkt efter avslutad mätning.





## 3.6 Programvara

#### 3.6.1 Introduktion/avsedd användning

Instrumentets programvara för kontroll och dataanalys, **Magellan**, levereras med instrumentet.

Magellan är en universell **programvara för styrning av läsare och dataanalys** för analysering av data som genereras av tester med mikroplattor med hjälp av mätinstrument från Tecan.

Magellan är tillgängligt i två versioner:

- Magellan och
- Magellan Tracker

Programvaran **Magellan** är avsedd för slutpunktsanalyser, kinetiska analyser och analyser med multipla inmärkningar med INFINITE F50 PLUS-instrument i enlighet med den avsedda användningen. Se kapitel 2.1 Avsedd användning.

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



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.

## 3.6.2 Installationprocedur

För att installera programvaran, sätt in USB-minnet i USB-porten och fortsätt på följande sätt:

- Magellans installationsguide startar automatiskt och guidar dig genom installationsprocessen. Om den inte gör det, kör filen "E:\Tecan.exe" (där E är USB-minnets enhetsbeteckning).
- 2. Välj "Magellan software" och beroende på vilken version du har beställt väljer du "Install Magellan" eller "Install Magellan Tracker" för att starta installationsproceduren och följ guiden.
- 3. Klicka på **Install** (Installera) för att starta installationsproceduren för programvaran.
- 4. Klicka på **I accept the terms of the license agreement** (Jag godkänner villkoren i licensavtalet) och **Next** (Nästa) för att fortsätta.
- 5. Sidan **Customer Information** (Kundinformation) visas: ange användarnamn och organisation.
- 6. Sidan Configuration (Konfiguration) visas: välj språk.
- 7. Sidan **Use for regulated environments** (Används i reglerade miljöer): klicka på **Next** (Nästa) för att fortsätta.
- 8. Klicka på Install (Installera) för att starta installationen.
- 9. Klicka på **Finish** (Slutför) för att avsluta installationen och stänga installationsprogrammet.

#### 3. Komma igång



Programvaran kan startas via Windows **Start**-menyn genom att välja **Magellan** i programgruppen Tecan.



i

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

Anmärkning Magellan V7.5 kan inte installeras tillsammans med andra versioner av Magellan-programvaran.

Som standard lagras alla filtyper som är associerade med Magellan i motsvarande underkataloger i följande katalog:

- Windows XP: C:\Documents och Settings\All Users\Documents\Tecan\Magellan
- Windows 7, Windows 8, Windows 10: C:\Users\Public\Documents\Tecan\Magellan

## 3.6.3 Installationskvalificering

Kontrollera att installationen av Magellan har slutförts med den automatiska programmet för installationskvalificeringen:

Starta *TecanIQ.exe* från standardinställningsvägen (C:\*Program Files\TecanWagellan)* eller från **Windows Start-menyn. Start > Programs > Tecan > MagellanIQ.** 

Klicka på **Check** (Kontrollera) för att börja installationskvalificeringen. Alla installerade komponenter bör ha status **OK**. Kontakta din lokala säljare för Tecan om eventuella problem rapporteras.

För att stänga installationskvalificeringsprogrammet, klicka på **Cancel** (Avbryt) eller **Exit** (Avsluta).



Anmärkning Installationskvalificeringen ska upprepas varje gång Magellan installeras eller uppdateras till en nyare version.



#### 3.6.4 Börja arbeta med Magellan

Den huvudsakliga typen av användargränssnittet i Magellan är guiden.

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 inmatningar kan göras.

#### Användargränssnitt – Guide-lista

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

Wizard List - Magellan		
$\frown$	What do you want to do?	
Start measurement	The Start Measurement wizard helps you to perform a measurement.	
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	•	
Exit Magellan		-

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 att 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.

Efter slutförd mätning skapas en arbetsyta.

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

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

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

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

#### 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.

#### 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.



Anmärkning För mer information om programvaran se Magellans användarhandbok.

Observera att vissa funktioner som beskrivs i Magellans användarhandbok kanske inte är relevanta (tillgängliga) i Magellan V7.5 i kombination med INFINITE F50 PLUS. All nödvändig information finns dock beskriven i denna användarhandbok.



Anmärkning Se detaljerat exempel på en ELISA-mätning i kapitel 6.

Sunrise fördefinierade metodfiler som skapats med Magellan V7.x eller senare kan öppnas med Magellan V7.5, mätparametrarna konverteras automatiskt.

Sunrise-metoder som skapats med tidigare versioner måste konverteras med alternativet "Convert From" (Konvertera från) i menyn övrigt/filhantering.



# 3.7 Magellan - Measurement Parameter Editor (Redigerare för mätparametrar)

**Measurement Parameter Editor** (Redigeraren för mätparametrar)används för att skapa arbetsflöden. Varje arbetsflöde skapas enkelt genom att dra och släppa processtegen i en sekvens enligt tillämpningen. Tillämpningens arbetsflöde är då synligt för användaren i panelen för arbetsflödet. Varje processteg (programelement) kan kopieras och klistras in (med hjälp av Windows standardgenvägar **Ctrl-C**, **Ctrl-V** eller den sammanhangsberoende musmenyn) och flyttas till den önskade positionen i arbetsflödet.

Measurement parameters - infinite F5			
🔶 Lab Ware 🔹	🔷 ▼ Plate	1	Selection
<ul> <li>Plate</li> <li>Part of Plate</li> </ul>	Plate definition: [GRE96t] - Greiner 96 Flat Transparent	Detais	Norning selected
Para a		Use a part of the plate	
Absorbance	Sector Plate A1:H12	2	
	🖕 🔻 Absorbance	3	
Accions	Wavelength Measurement: 405 mm		
Move Plate	Reference: 405 rm		
😡 Kinetic 🔹			
C Kinetic Cycle Kinetic Condition			
Miscellaneous 🛞			
E Comment			
Wait (Timer)			
Jincubation			
Kontrollfölt	Donal för arbeteflöda		
Kontrolliali	Faller for arbeishoue		Info
			pane
		$\sim$	•
		202	
			×
Number of plates: 1			
Help		OOCE MEACUDEMENT D. C.	
Cancel <<< Back	СН	UUSE MEASUREMENT PARAMET	ERS 🕑
© 2009 Tecan			

**Measurement Parameter Editor** (Redigeraren för mätparametrar) består av följande alternativ som beskrivs i detalj i följande kapitel:

- Kontrollfält
- Panel för arbetsflöde
- Infopanel



#### 3.7.1 Kontrollfält

Kontrollfältet är indelad i fem sektioner. Varje avsnitt innehåller programelement som används för att skapa ett individuellt arbetsflöde.

Skapa ett arbetsflöde antingen genom att dubbelklicka på det valda programelementet eller genom att dra och släppa det i panelen för arbetsflöde.

Följande programelement är tillgängliga när du använder INFINITE F50 PLUS:

Lab Ware (Laboratoriearti klar)	Plate (Platta) Part of Plate (Del av platta)
Measurements (Mätningar)	Absorbance (Absorbans)
Actions (Åtgärder)	Shaking (Skakning) Move Plate (Flytta platta)
Kinetic (Kinetik)	Kinetic Cycle (Kinetikcykel) Kinetic Condition (Kinetikförhållande)
Miscellaneous (Övrigt)	Comment (Kommentar) User Request (Användarförfrågan) Wait (Timer) (Vänta - timer) Incubation (Inkubation)

#### Lab Ware (Laboratorieartiklar)

#### Plate (Platta)

Programelementet **Plate** (Platta) används för att välja ett format för plattan från rullgardinsmenyn **Plate definition** (Plattdefinition). Klicka **Details**... (Detaljer) för att se ytterligare information om den valda plattan.

🔷 🔻 Plate		1
Plate definition:	[COS96it] - Corning 96 Flat Transparent	Details
		Use a part of the plate

#### Part of Plate (Del av platta)

Som standard är programelementet **Part of Plate** (Del av platta) infällt. När den utvidgas genom att klicka på visas en mikroplatta med 96 brunnar. För att mäta enskilda brunnar klickar du på den önskade brunnen eller om du vill mäta ett intervall av brunnar markerar du det önskade området. När du klickar på **Details...** (Detaljer), kan du zooma in på förhandsgranskningen av plattan.

🔹 🔻 Part of Plate		2
1       2       3       4       5       6       7       8       9       10       11       12         B       0 <th>Details</th> <th></th>	Details	



#### Individuella delar av platta

Individuella delar av plattan kan väljas:



Ett andra intervall av brunnar kan väljas genom att trycka på **Kontrolltangenten** på tangentbordet och markera brunnar som ska väljas.

#### Measurements (Mätningar)

#### Absorbance (Absorbans)

Programelement **Absorbance** (Absorbans) används för att utföra absorbansmätningar. Ange eller välj respektive parametrar.

Två rullgardinsmenyer visar tillgängliga våglängder för mät- och referensfilter enligt de insatta absorbansfiltren. Om rullgardinsmenyerna är tomma har filtren inte definierats.

🍦 🔻 Absorb	ance			3
Wavelength-	400		Label	
Measurement:	492 nm	*	Name: Labell	
Reference:	405 nm	~		

#### Actions (Åtgärder)

#### Shaking (Skakning)

Välj programelementet **Shaking** (Skakning) om plattan ska skakas, antingen före mätningen eller mellan kinetikcykler.

📉 🔻 Shaking		4
Parameter Duration: 1 📚 sec	Intensity: Wide Amplitude: 14,1 mm Frequency: 2,1 Hz	Wait a couple of seconds

Ange respektive parametrar.

Duration (Varaktighet)	Ange skakningens varaktighet.
Intensity (Intensitet)	Ange önskat skakningsläge. Amplitud och frekvens visas när du väljer respektive skakningsläge.

Se kapitel 4.1.1 Skakning av mikroplatta för tillgängliga skakningslägen.



Om du klickar på länken <u>Wait a couple of seconds</u> (Vänta ett par sekunder) läggs ett nytt programelement till. Se sida 25 för mer information.

**Kinetic (Kinetik)** 

#### Kinetic Cycle (Kinetikcykel)

Använd programelementet **Kinetic Cycle** (Kinetikcykel) för att utföra flera följande mätningar som kan utföras inom vissa intervaller.

🎧 🔻 Kinetic Cycle		3
Ovcles       Image: Second state       Duration	Kinetic Interval         Image: Use kinetic interval.         Time:       00:01:00 Image: [hh::mm:ss]         Time:       60000 Image: ms	

Ange respektive parametrar.

Cycles (Cykler)	<ul> <li>Number of cycles (Antal cykler): Ange ett nummer eller klicka på pilarna uppåt eller nedåt för antalet faktiska mätsteg (2–1000 cykler).</li> <li>Duration (Varaktighet): Ange varaktighet, format hh:mm:ss.</li> </ul>
Kinetic Interval (Kinetikintervall)	<b>Use kinetic interval</b> (Använd kinetikintervall): Ange tidsintervall (hh:mm:ss eller ms).

#### Kinetic Condition (Kinetikförhållande)

Använd programelementet **Kinetic Condition** (Kinetikförhållande) för att definiera vilka åtgärder som ska utföras i en viss cykel.



Om **2** anges för **Execute command at cycle** (Utför kommandot vid cykeln) inom en kinetisk mätning som kan innehålla steget **Shake** (Skakning), utförs endast skakning i cykel 2.



Anmärkning Kinetikhållanden som Shake (Skakning) bör sättas in direkt efter programelementet Kinetic Cycle (Kinetikcykel) för att säkerställa optimal reproducerbarhet av resultaten. Användarna rekommenderas att ställa in lämpliga skript innan mätningarna och att använda samma skript för alla liknande kinetiska

mätningar för att få jämförbara resultat.



#### Miscellaneous (Övrigt)

#### **Comment (Kommentar)**

Använd programelementet **Comment (Kommentar)** för att skriva in en anmärkning eller ett uttalande för den aktuella mätningen i textfältet.

🚛 🔻 Comment	6
Comment:	

#### User Request (Användarförfrågan)

Programelement **User Request** (Användarförfrågan) informerar instruments operatör att utföra en bestämd åtgärd under arbetsflödet vid en viss tidpunkt.

🚯 🔻 User Request	7
Text:	

Om exempelvis programelementet **Move Plate** (Flytta platta) används för att flytta plattan för att utföra en viss åtgärd, ska den inskrivna texten informera operatören om att utföra dessa åtgärder. En dialogruta visar meddelandet och mätningen stannar tills du klickar på **OK**.

#### Wait (Timer) (Vänta - timer)

Använd programelementet **Wait (Timer)** (Vänta - timer) för att definiera en viss väntetid innan nästa steg i ett arbetsflöde utförs.

Ange önskad tid i fältet Wait time (Väntetid).

🥝 🔻 Wait (Timer)		5
Timer Wait time: 00:01:00 ♥ (hh::mm:ss)	Options Wait for injection Ignore wait at last kinetic cycle	

Ange respektive parametrar:

Timer	Ange Väntetiden (hh:mm:ss)
<b>Options</b> (tillval)	Ignore wait at last kinetic cycle: (Ignorera väntan vid den sista kinetikcykeln:) När programsteget Vänta (Timer) är den sista åtgärden i en kinetisk omgång kommer väntetiden att ignoreras i den sista cykeln.

#### **Incubation (Inkubation)**

🥹 ▼ Incubation	4
Timer Incubation time: 00:01:00 💸 (hh:mm:ss)	

Ange lämpliga parametrar för inkubation:

Incubation time	Ange den totala tiden (min. 5 s)
(Inkubationstid)	



## 3.7.2 Workflow Pane (Panel för arbetsflöde)

**Panelen för arbetsflöde** är fönstret där mätskriptet är synligt och parametrar definieras och redigeras.

Det finns två sätt att lägga till ett programelement från Kontrollfältet till **Panelen** för arbetsflöde:

- Välj ett programelement från Kontrollfältet, genom att dubbelklicka på det läggs det till i Panelen för arbetsflöde direkt efter det föregående programelementet.
- Klicka på programelementet i Kontrollfältet och dra det till Panelen för arbetsflöde till respektive position.

Programelementen är numrerade enligt deras ordningsföljd.

När ett programelement har lagts in i **Panelen för arbetsflöde** kan inställningar och parametrar för detta element anges eller redigeras.

Enstaka programelement i **Panelen för arbetsflöde** kan fällas ihop för att visa den viktigaste informationen eller expanderas för att få tillgång till alla funktioner som kan redigeras. Klicka på en av trianglarna bredvid programelementets titel, eller , för att växla mellan de två visningslägena.

Som standard börjar redigeraren av mätparametrar med elementet **Plate** (Platta), elementet **Part of Plate** (Del av platta) (hopfällt) och ett **Absorbance** (Absorbans)-element i **Panelen för arbetsflöde**.

De aktuellt valda programelementen i **Panelen för arbetsflöde** visas med en gul linje i den övre kanten.

Om ett programelement innehåller fel eller är ogiltigt i det aktuella arbetsflödet flaggas elementet med en felmarkering och elementets nummer markeras med rött. I **Infopanelen** visas detaljerad information om felet. Om arbetsflödet innehåller fel kan mätparametrarna inte väljas.

#### Elementens hierarki

Elementens hierarki i Panelen för arbetsflöde är följande:

- 1. Plate (Platta)
- 2. Part of Plate (Range) (Del av platta intervall)

Alla önskade mätsteg kan läggas till direkt efter ett platt- eller intervall-element. Använd **Release** (Släpp) och **Indent** (Dra in) för att ändra sekvensen för utförandet av den enskilda stripskomponenten. Välj ett element i **Panelen för arbetsflöde**, klicka på höger musknapp och välj **Release** (Släpp) eller **Indent** (Dra in).

Andra element från **Kontrollfältet** kan läggas in i ett arbetsflödes hierarki på följande sätt:

Det första **Range** (Intervall)-elementet läggs till direkt efter elementet **Plate** (Platta), därefter kan alla efterföljande **Range** (Intervall)-element läggas till.

Kinetiska steg är möjliga inom elementen Plate (Platta) eller Range (Intervall).

Stegen **User Request** (Användarförfrågan), **Comment** (Kommentar) och **Wait** (Vänta) är möjliga inom elementen Plate (Platta) eller **Range** (Intervall).

#### 3.7.3 Infopanel

**Infopanelen** på höger sida av skärmen visar information som är relevant för det markerade programelementet. Eventuella varningar och fel visas.



# 3.8 Magellan - Definiera mätningar

I följande kapitel beskrivs några exempel för att illustrera definitionen av olika mätningar.

#### 3.8.1 Definiera slutpunktsmätningar

Följande exempel beskriver en **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.

🔷 🔻 Plate		1
Plate definition: [COS96/t] -	Corning 96 Flat Transparent	Use a part of the plate
	6 7 8 9 10 11 12 Details	2
↓ ▼ Absorb Wavelength Measurement	450 mm	Label1



## 3.8.2 Definiera mätningar med multipla inmärkningar

Följande exempel beskriver en **Mätning med multipla inmärkningar för absorbans** i ett definierat intervall i en mikroplatta med 96 brunnar (A1:E7). 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. Klicka på för att expandera elementet **Part of Plate** (Del av platta). Välj sedan det önskade plattintervallet (A1:E7).
- 3. Ange önskad mätningsvåglängd.
- 4. Lägg till 2 eller fler **Absorbance** (Absorbans)-element och ange mätningsvåglängderna.

🔷 🔻 Plate						1
Plate definition:	[COS96ft] - Corning 96 Flat Transparent				~	Details
					<u>Use a p</u>	art of the plate
🄹 🔻	Part of Plate					2
A B C D E G H		Details				
4	▼ Absorbance					3
٢٧	Vavelength	Label		_		
M	feasurement: 405 nm 👻	Name:	Label1	•		
	Reference: 405 nm					
4	▼ Absorbance					4
-V	Vavelength	Label				
M	feasurement: 492 nm 💙	Name:	Label2	<b>v</b>		
	Reference: 405 nm					
4	▼ Absorbance					5
- ۷	Vavelength	Label				
M	leasurement: 620 nm 👻	Name:	Label3	•		
	Reference: 405 nm					
						63



#### 3.8.3 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. Lägg till 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.

🔶 🔻 Plate	,	1
Plate definition:	[COS96/t] - Corning 96 Flat Transparent	Details Use a part of the plate
	Part of Plate	2
1 2 A 0 C 0 C 0 C 0 C 0 C 0 C 0 C 0 C	2 3 4 5 6 7 8 9 10 11 12	
Ģ	🖓 🔻 Kinetic Cycle	3
	Cycles       Similar Cycles:       50 ♀         O Duration       Image: Cycles:       100:02:30 ♀         (http://www.ss)       Time:       150000 ♀         Time:       150000 ♀       ms	
	🐇 🔻 Absorbance	4
	Wavelength Measurement: 492 nm V Reference: 405 nm V	



#### 3.8.4 Dra in och släppa programelement

Beslutet att dra in/släppa ett programelement kommer att ändra instrumentets arbetsflöde under mätningar.

Åtgärderna i alla programelement med samma indragning utförs i sekvens. Det enda sambandet mellan dessa programelement är att nästa åtgärd börjar direkt efter att den föregående åtgärden är slutförd.

Ett programelement som är mer indraget än det föregående programelementet visar att de två programelementen är beroende av varandra. Detta innebär att de parametrar som definieras i det första programelementet också är aktiva för det andra (indragna) programelementet.

Följande är ett exempel på hur man definierar en **Kinetik med multipla** inmärkningar med två **Absorbansmärkningar**. Exemplet visar att de två programelementen **Absorbance** (Absorbans) är beroende av programelementet **Kinetic Cycle** (Kinetikcykel), som är beroende av programelementet **Part of Plate** (Del av platta), som är beroende av programelementet **Plate** (Platta). Definiera parametrarna för ett exempel på följande sätt:

- 1. Plate (Platta): e.g., Greiner 96 Flat Transparent
- 2. Kinetic Cycle (Kinetikcykel)/Number of cycles (Antal cykler): 5
- 3. Absorbance/Wavelength Label 1 (Absorbans/våglängdsmärkning 1): 450 nm
- 4. Absorbance/Wavelength Label 2 (Absorbans/våglängdsmärkning 2): 492 nm

Panelen för arbetsflöde visas enligt skärmdumpen:

🔷 🔻 Plate			1
Plate definition: [GRE96ft] - 0	åreiner 96 Flat Transparent		Use a part of the plate
🔹 🔻 Part of Plate			2
1       2       3       4       5       6         B       ••••••••••••••••••••••••••••••••••••	7       8       9       10       11       12         0 <th>ils</th> <th></th>	ils	
😡 🔻 Kinetic (	Cycle		3
© Number of cy O Duration	cles: 5	⊂ Kinetic Interval	
🕴 🕶 At	sorbance		4
Waveleng Measurem Referen	th ent: 450 nm 💌 nce: 405 nm 😪	Label Name: Label1	
🖕 🔫 At	sorbance		5
Waveleng Measurem Referen	th ent: 492 nm 💌 ice: 405 nm 😪	Label Name: Label2	
			10 m

Definitionen ovan resulterar i följande arbetsflöde:

Absorbansen av alla brunnar i mikroplattan med 96 brunnar mäts först vid 450 nm och sedan vid 492 nm. Båda absorbansmätningar utförs i fem kinetikcykler.



Genom att släppa det andra programelementet **Absorbance** (Absorbans) så att det är i linje med elementet **Kinetic Cycle** (Kinetikcykel) förändras arbetsflödet. Välj det andra programelementet **Absorbance** (Absorbans) och klicka på höger musknapp. Välj **Release Strips** (Släpp strips) från den sammanhangsberoende menyn. **Panelen för arbetsflöde** visas enligt följande skärmdump:

Plate definition:	[GRE96lt] - Greiner 96 Flat Transparent		Use a part of the plate
🤹 🔻 Pa	rt of Plate		2
1 2 A • • • • C • • • D • • • F • • • G • • H • • •		)etais	
	▼ Kinetic Cycle		3
©	les Number of cycles: 5	Kinetic Interval	
	💡 🔻 Absorbance		4
	Wavelength Measurement: 450 nm	Label Name: Label1	
	▼ Absorbance		5
Wa Mea	velength asurement: 492 nm 💌 Reference: 405 nm 👻	Name: Label	
			K

l detta arbetsflöde görs en **Absorbanskinetisk** mätning med 5 cykler för den första absorbansen vid 450 nm, efter denna utförs **Slutpunktsmätning för absorbans** vid 492 nm.



## 3.9 Optimering för bästa prestanda

Instrumentet har fabrikstestats för att säkerställa att dess prestanda ligger inom de angivna gränserna (se 4.4.2 Mätspecifikationer för detaljer).

Den största noggrannheten kan erhållas från instrumentet genom att observera de rekommendationer som nämns nedan.

#### 3.9.1 Instrumentplats

Instrumentet ska placeras på en lämplig plats (se kapitel 3.3 Miljökrav för detaljerad information).

#### 3.9.2 Tillvägagångssätt

#### Allmänt

- Det rekommenderas att man följer de standardiserade tillvägagångssätt för de analyser som används.
- Den bästa reproducerbarheten uppnås när mätningsvåglängden motsvarar den maximala absorbansvåglängden för den särskilda lösningen. Det är viktigt att använda den maximala absorbansvåglängden om provets absorbanskurva ligger på ett smalt våglängdsband. Observera att mätningar av lutningen på en absorbanstopp begränsar noggrannheten hos OD-värdena.
- Efter mätning av varje mikroplatta, se provsatsens förpackning för information om valideringsproceduren
- Använd de rekommenderade absorbansfiltren för INFINITE F50 PLUS.

#### **Mikroplattor**

 Instrumentet kan användas med de typer av mikroplattor som beskrivs i kapitel 4.4.3 Mikroplattor. De bästa resultaten uppnås när man använder mikroplattor med plan botten. Mätresultaten kan variera beroende på vilken typ av mikroplatta som används.
 Var särskilt försiktig när du använder mikroplattor med C-, U- eller V-

formad botten eller strips eftersom det är möjligt att mätresultaten kan skilja sig från de specifikationer som beskrivs i detta dokument. Se till att typen av mikroplatta som används med INFINITE F50 PLUS absorbansläsare är lämplig för respektive tillämpning.

- Använd endast helt rena mikroplattor.
- Låt inte damm lägga sig på lösningarna eller mikroplattan under inkubationsperioden innan mätningen. Det rekommenderas att använda ett skydd när en mikroplatta placeras utanför instrumentet.
- Felaktigheter i mängden lösning som pipetteras har en större effekt på resultat som erhålls när små mängder lösningar används.
- Formen på lösningens menisk kan orsaka felaktigheter i resultaten, särskilt om små mängder lösning används.



# 4. Instrumentfunktioner

# 4.1 Instrumentfunktioner

Följande lägen för absorbansmätning är tillgängliga på INFINITE F50 PLUS:

slutpunktsmätningar, kinetiska mätningar och mätningar med multipla inmärkningar.

#### 4.1.1 Skakning av mikroplatta

The INFINITE F50 PLUS kan skaka mikroplattan innan mätning. Mikroplattan kan också skakas mellan varje kinetiska mätcykel.

Använd Magellan för att ställa in skakningsläget.



#### FÖRSIKTIGHET NÄR MIKROPLATTOR SKAKAS KAN SPILL UPPSTÅ OM BRUNNARNA ÖVERFYLLS.

Skakningsläge för INFINITE F50 PLUS:

Skakläge	Skakbredd	Skakfrekvens
HIGH (HÖG)	2,8 mm	12,3 Hz
NORMAL	4,4 mm	9,2 Hz
LOW (LÅG)	4,4 mm	7,8 Hz
BRED	14,1 mm	2,0 Hz

# 4.2 Instrumentbeskrivning

Illustrationen nedan visar instrumentets komponenter.





#### 4. Instrumentfunktioner

Status-LED ger information om instrumentets status:

- Grönt blinkande: instrumentet är inte anslutet till Magellan
- Grön: instrumentet är anslutet och redo för mätning
- Röd: mätning pågår

På instrumentets vänstra panel finns USB-porten, huvudströmbrytaren och eluttag.



Typskylten finns på instrumentets undersida.

Exempel på typskylt



Innehållet på typskylten (t.ex modellnamn och artikelnummer) kan variera beroende på specifik modell.

Se "Declaration of Conformity" på sista sidan i detta dokument för en översikt över de olika instrument som denna användarhandbok gäller för.

# 4.3 Beskrivning av filterhjul

INFINITE F50 PLUS standardfilterhjul levereras med fyra smalbandsinterferensfilter som har en fast våglängd (405, 450, 620 och 492 nm). Det är möjligt att utrusta filterhjulet med upp till åtta filter. Kontakta din lokala säljare för Tecan för tillgängliga filter.

Filtren i standardfilterhjulet monteras enligt nedanstående:

Filterposition	Filtervåglängd
1	405 nm
2	450 nm
3	620 nm
4	492 nm
5–8	tomma filterposition



När en våglängd väljs för mätning förs det specifika filtret in i ljusstrålen genom att flytta filterhjulet till rätt position.



Anmärkning För mer information om definition av ett nytt filter, se 7.5.2 Definiera filter.



# 4.4 Instrumentspecifikationer

I tabellen nedan anges specifikationerna för absorbansläsaren INFINITE F50 PLUS.

# 4.4.1 Allmänna specifikationer

PARAMETRAR	EGENSKAPER		
Strömingång Extern strömförsörjning	Strömförsörjning Grundläggande instrumentet med AC-adapter: 100 - 240 V AC, 50/60 Hz, max. 1.2 A (automatisk avkänning, överspänningskategori II) Grundläggande instrumentet utan AC-adapter: 24 V DC (överspänningskategori I)		
Förbrukning INFINITE F50 PLUS	Standby-läge: ca. 12 VA Driftläge: max. 30 VA		
Yttermått	Bredd: 34,7 cm (13,66 inch) Djup: 18,9 cm (7,44 inch) Höjd: 13,4 cm (5,28 inch)		
Vikt	2,6 kg (strömförsörjning inkluderad)		
Omgivningstemperatur:			
Drift	15 °C till 35 °C (59 °F till 95 °F)		
Förvaring	-30 °C till 60 °C (-22 °F till 140 °F)		
Relativ luftfuktighet	20 % till 80 %		
Föroreningsgrad	2		
Avfallshantering	Kontaminerat avfall		
Miljö	Se kapitel 3.3 Miljökrav för mer information.		


# 4.4.2 Mätspecifikationer

PARAMETRAR	EGENSKAPER					
Mättid: enkel våglängd dubbel våglängd	< 15 sekunder < 20 sekunder					
Våglängdsintervall: Standard	400–750 nm					
Mätningsintervall: 400–750 nm	0–4,000 OD					
Upplösning:	0,0001 OD					
Noggrannhet: 450, 492 nm 0,000–2,000 OD 2,000–3,000 OD	≤ (1,0 % + 0,010 OD)* ≤ (1,5 % + 0,010 OD)*					
Precision: 450, 492 nm 0,000–2,000 OD 2,000–3,000 OD	≤ (0,5 % + 0,005 OD)* ≤ (1,0 % + 0,005 OD)*					
Linjäritet: 450, 492 nm 0,000–2,000 OD 2,000–3,000 OD	≤ 1 % ≤ 1,5 %					
Våglängdsval: Standardfilter	Smalbandsinterferensfilter. Upp till åtta filter kan monteras i ett filterhjul.					
Noggrannhet filtervåglängd:	Central våglängd ± 2 nm					
Filterbandbredd: Vid 50 % transmission	10 ± 2 nm					
Ljuskälla:	LED					
Datorgränssnitt:	USB					
All ansluten utrustning måste vara godkänd enligt IEC 60950-1 – Utrustning för						

informationsbehandling – Säkerhet och motsvarande lokala standarder.

\* bättre än eller lika med x % av mätvärdet plus tillhörande OD-värde.

## 4.4.3 Mikroplattor

Alla mikroplattor med 96 brunnar med transparent botten (plan, C-, U- och Vformade, inklusive strips) som överensstämmer med följande standarder kan användas med INFINITE F50 PLUS absorbansläsare:

ANSI/SBS 1-2004; ANSI/SBS 2-2004; ANSI/SBS 3-2004; ANSI/SBS 4-2004



#### FÖRSIKTIGHET ANVÄND ENDAST MIKROPLATTOR UTAN LOCK OCH INTE MIKROPLATTOR SOM ÄR HÖGRE ÄN 15,2 MM.



## Hantering av mikroplattan

Sätt in eller ta ut mikroplattan endast när platthållaren är helt utskjuten (enligt illustrationen nedan) och platthållarens motor inte är aktiv. Öppna inte höljets lock när LED-statusen lyser rött.



### VARNING BÄR ALLTID ENGÅNGSHANDSKAR OCH SKYDDSKLÄDER NÄR DU HANTERAR MIKROPLATTAN.



Placering av brunn A1 Platthållare – helt utskjutet

# 4.5 Instrumenttillbehör

Listan nedan innehåller de tillgängliga valfria tillbehören för INFINITE F50 PLUS, som kan beställas:

- Extra filter
- Verktyg för filtersammansättning
- MultiCheck<sup>TM</sup>-plattan för serien INFINITE F50

För mer information och tillgänglighet i ditt land, kontakta din lokala säljare för Tecan.



# 5. Kvalitetskontroll

# 5.1 Inledning



#### FÖRSIKTIGHET OM INFINITE F50 PLUS ANALYTISKA PRESTANDA VID NÅGOT TILLFÄLLE KAN IFRÅGASÄTTAS SKALL ANVÄNDAREN FÖLJA DE ANVÄNDARINSTRUKTIONER SOM GÄLLER FÖR KVALITETSKONTROLL ELLER KONTAKTA TECANS LOKALA KUNDSUPPORT.

Det här kapitlet tillhandahåller information om självkontroll för instrumentet och instruktioner om hur du enkelt kan kontrollera driftskvaliteten.

# 5.2 Självkontroll

Under anslutningen av INFINITE F50 PLUS till Magellan läsarens kontrollprogramvara kontrolleras motorer och sensorer, och platthållaren och filterhjulet initialiseras.

Före varje mätning utförs en självkontrollkalibrering för att säkerställa att instrumentet fungerar korrekt och för att kalibrera det optiska systemet.

# 5.3 Driftskvalificering (OQ)

Följande test kan göras för att säkerställa att instrumentet fungerar korrekt och att exakta resultat erhålls.

Instrumentets repeterbarhet och noggrannhet kan variera beroende på vilken typ av lösning och mikroplatta som används.

För att eliminera denna effekt fabrikstestas instrumenten med en kalibreringsplatta, som tar bort lösningens inverkan och eventuella variationer beroende på mikroplattans placering vid mätningen.

# 5.3.1 Tester med MultiCheck

Tester med MultiCheck tillhandahåller en automatiserad kontroll av läsarens prestanda inklusive noggrannhet, linjäritet, precision och inriktning med standarder som kan påvisas med NIST.

# 5.3.2 Tester med mikroplattor

Om de optiska densiteterna i mikroplattans brunnar inte är konsekventa kommer de resultat som erhålls med denna typ av mikroplatta att påverkas.

Denna inkonsekvens kan kontrolleras genom avläsning av en tom mikroplatta.

De OD-värden som erhålls vid mätningen av den tomma mikroplattan bör ligga inom ett smalt intervall. Till exempel:  $\pm 0,010$  OD.

Om OD-värdena inte ligger inom detta intervall bör denna typ av mikroplatta inte användas.

Genom användning av mätningar med dubbla våglängder tas skillnaderna i mikroplattans OD-värden bort eller reduceras till en nivå som ligger inom godtagbara gränser.



## 5.3.3 Instrumentprecision med flytande prover

Denna procedur kan användas för att kontrollera mätningarnas reproducerbarhet. Användning av en mikroplatta med plan botten rekommenderas.

Fyll en ny mikroplatta med nypreparerad Orange G-lösning. Använd olika spädningar av lösningen i varje brunn så att ett intervall av optiska densiteter erhålls. Se till att brunnarna innehåller minst 200 µl. Spädningsserien bör vara inom 0,1 till 3,0 OD. För att nå ca 3 OD rekommenderas 125 mg.l<sup>-1</sup>Orange G (Sigma, Cat. Nr. 07252).

Programmera ett test med 492 nm filter och mät sedan mikroplattan minst tre gånger.

Beräkna följande för varje brunn:

- genomsnittliga OD-värdet
- standardavvikelse

### Exempel

#### Avläsningar 0,000 till 2,000 OD

Standardavvikelsen för varje brunn bör ligga inom (0,5 % + 0,005 OD).

Beräkning av högsta tillåtna avvikelse med 1,000 OD som genomsnittligt ODvärde:

1,000 \* 0,5 % + 0,005 = 0,010 OD

#### Avläsningar 2,001 till 3,000 OD

Standardavvikelsen för varje brunn bör ligga inom (1,0 % + 0,005 OD).

Beräkning av högsta tillåtna avvikelse med 2,400 OD som genomsnittligt ODvärde:

2,400 \* 1,0 % + 0,005 = 0,029 OD

#### Avläsningar över 3,000 OD

Avläsningar över 3,000 OD används endast som en indikering och precisionen kan inte garanteras.

## 5.3.4 Instrumentlinjäritet med flytande prover

Instrumentets linjäritet och tillämpning vid den använda våglängden kan kontrolleras med hjälp av en spädningsserie för en lösning.

Resultatet beror på renheten hos den använda färgen och menisken hos vätskan i brunnarna.

Som referens kan en spädningsserie av Orange G-lösning för mätningar vid 492 nm användas.

Spädningsserien bör vara inom 0,1 till 3,0 OD. För att nå ca 3 OD rekommenderas 125 mg.l<sup>-1</sup>Orange G (Sigma, Cat. Nr. O7252).

För andra våglängder måste olika lösningar användas.

En volym av 200 µl av varje spädning pipetteras på mikroplattan. Ett minimum av minst två prover bör användas för varje spädning för att reducera de fel som orsakas av pipetteringen.



Mikroplattan mäts sedan och en linjär regression av OD mot koncentration ritas med utgångspunkt från medelvärdet av de uppmätta OD-värdena.

Beräkna det oviktade residualkvadratsumman R<sup>2</sup> från regressionslinjen.

Typiska residualkvadratsummor för en standardapplikation är lika med eller bättre än  $R^2 = 0,998$ .



Anmärkning Data kan variera på grund av felaktigheter vid pipetteringen.



# 6. Exempel på tillämpning

# 6.1 Inledning

Magellans **exempelfiler** tillhandahåller Magellan-metoder och arbetsytor som introducerar programmet och underlättar användarens arbete med det. Exempelfilerna för en kvantitativ och en kvalitativ ELISA-analys installeras automatiskt vid installation av Magellan.

# 6.2 Steg-för-steg exempel: Kvantitativ ELISA

I det här kapitlet finns ett steg-för-steg exempel (kvantitativt test) på hur man skapar en metod i Magellan. Genom att följa instruktionerna lär du dig att definiera utvärderingar från en provsatsbeskrivning i Magellan.



Anmärkning Exempelfiler visas automatiskt i metodlistan i Magellan. Dessa filer finns tillgängliga för Magellan Tracker i standardsökvägen för data och måste konverteras.

# 6.2.1 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									
E	C1	C5	S3									
F	C2	C6										
G	C2	C6										
Н	C3	S1										
	05	51										

BLK = Blank, NC = Negativ kontroll, C1 - C6 = Kalibratorer (Standarder), S1 - S... = Prover



### Mätning och utvärdering

Läs av plattan vid våglängden 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–OD 620) 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 (arbetsyta) automatiskt och en rapport med mätparametrar, plattlayout, blanka värden, standardkurva, IgM-koncentrationer, definition av cutoff, kvalitativa resultat av proverna och valideringar skapas.

Dessutom lagras layouten och de kvalitativa resultaten som ASCII-filer.



# 6.2.2 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 **Continue** (Fortsätt) på **Välkomstsidan** av **Create/edit a method wizard** (Guiden Skapa/redigera en metod) och dialogrutan **Select a file** (Välj en fil) visas. Välj **New** (Ny).

Create/Edit a Method			
O New ⊙ Open Commth	Show:	Files from this instrument	Print Preview Status
		_	
Help Cancel <<< Back		MAKE YOUR S	
		© 2	009 Tecan

#### Mätparameterar

Klicka på **Make Your Selection** (Gör ditt val) och sidan **Measurement parameters** (Mätparameterar) visas:

Measurement parameters - infinite F5	0		×
🔷 Lab Ware 🔹	- Velate	1 Selection	
🔶 Plate	Plate definition: [GRE96t] - Greiner 96 Flat Transparent	Details     Nothing selected	
Part of Plate		Use a part of the plate	
Measurements (*)	Sector A1:H12	2	
4 Absorbance	Abardance		
Actions	Wavelenth		
🔀 Shaking	Measurement: 405 nm 💙 Name: Label1 🗸		
🐔 Move Plate	Reference: 405 nm		
Kinetic (*			
G Kinetic Cycle			
🏂 Kinetic Condition			
Miscellaneous 🛞			
E Comment			
<ul> <li>User Request</li> <li>Wait (Timer)</li> </ul>			
Incubation			
		203	
			~
Number of plates: 1	,	,	
Heb			
Cancel KKK Back	СНО	DOSE MEASUREMENT PARAMETERS	
			,



#### 6. Exempel på tillämpning

På **Wavelength** (Våglängd)-stripsen välj 492 nm som mätningsvåglängd och 620 nm as referensvåglängd.

🖌 🔻 Absorbance			3
Wavelength Measurement: 492 nm	V Label Name:	Label1 💌	

Fortsätt med guiden genom att klicka på **Choose measurement parameters** (Välj mätparametrar) och fönstret för **Plate layout (Plattlayouten)** visas.

Create/Edit a Method	10%	100% Zo	om											
Method layout														
	0	1	2	3	4	5	6	7	8	9	10	11	12	
												Identifiers	Exp. group:	
	B											ST PC NC	ID-Num	
	C											LPC HPC BF	Replicates	
	D											BF Def. identi		
	E											Fill selection	Delete selection	
	F													
	G													AN Austria
	H													TEC
Help Cancel <<< Back													NEXT D	



## **Utforma layout**

Definiera plattlayouten med hjälp av dialogrutan **Well Assignment** (Tilldelning av brunn) på höger sida av skärmen.

We	Well Assignm.								
r.	Identifier	s	Exp. group:						
	SM	^	1						
	BL		- ID-Num.						
	ST		≥ 4						
	PC								
	NC								
	LPC		- Replicates -						
	HPC		≠ 4						
	BF		• Fix number:						
	RF	~	1						
l	Def. ide	entif.	O All						
	Fill selecti	on	Delete selection						

I grupprutan Identifiers (Identifierare) välj BL (Blank).

I rutan Experimental group (Försöksgrupp) kvarstår numret 1.

I grupprutan Replicates (Replikat), väljs All (Alla) automatiskt.

Klicka på brunn A1, som då markeras med en röd kant.

Klicka på **Fill selection** (Fyll urval) och brunnen märks med den utvalda identifieringstypen.



#### Anmärkning En enskild 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)

l rutan Experimental group (Försöksgrupp) kvarstår numret 1.

All (Alla) i grupprutan Replicates (Replikat), väljs automatiskt.

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.

Kalibratorer (standarder) måste tilldelas till brunnar **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 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 valda brunnarna till som replikat till de befintliga replikaten. Med alla andra identifieringstyper läggs de valda 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.

l 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.

I 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 alternativet **Transformed data** (Transformerad data), för att definiera blankreducering.

En dialogruta kommer att fråga dig om du vill definiera en blankreducering. Klicka på **Yes** (Ja). Följande fönster visas:

Crea	te/Edit a Method															×
Un	do Redo		10%	100% Za	om											
	(	Ś	Input data	Difference data				~							Constants	Options
	Method layout	ę		fx xBL1				Line			~	Available dat	a (multiple data	sets) 🔻	Functions&Co	nstants 🔻
8	Plate layout	eate,														
	Conc, Dil, Refvalues	edit														
-	Transformed data 🛛 🔗	meth			-	-			-	_	_	_				$\sim$ 1
	Blank reduction	ā		) 1	2	3	4	5	6	7	8	9	10	- 11		
	Add new transformation	r			1	-					_	~			_	- 11
	Concentrations 🛞		0	BL1	ST1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37	- 11
	standard curve			x-BI 1	2/2 x-BI 1	Z/Z x-BL1	x-BL1	2/2 x-BL1	2/2 x-BL1	2/2 x-BL1	2/2 x-BL1	2/2 x-BL1	2/2 x-BL1	2/2 x-BL1	ZIZ x-BL1	- 11
-0	Evaluate data 🛞			NC1	ST1 4	SM1 2	SM1 6	SM1 10	SM1_14	SM1 18	SM1 22	SM1 26	SM1 30	SM1 34	SM1 38	- 11
	QC Validation			1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	- 11
100	Data has the o			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	- 11
	Data export			NC1	ST1_4	SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	- 11
8	Printed report			<u> </u>	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	- 11
8	Automated data handling			X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	- 11
8	Miscellaneous		ll r	SI1_1 1/2	SI1_5 1/2	SM1_3	SM1_/	SM1_11 1/2	SM1_15 1/2	SM1_19 1/2	SM1_23	SM1_27	SM1_31 1/2	SM1_35	SM1_39	- 11
	User prompts Number format			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	- 11
	Method notes			ST1 1	ST1 5	SM1 3	SM1 7	SM1 11	SM1 15	SM1 19	SM1 23	SM1 27	SM1 31	SM1 35	SM1 39	- 11
				2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	- 11
				x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	- 11
				ST1_2	ST1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	- 11
			J	=> 1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	
				ST1 2	ST1 6	SM1 4	SM1 8	SM1 12	SM1 16	SM1 20	SM1 24	SM1 28	SM1 32	SM1 36	SM1 40	.e
				2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	Aust
				x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	AND
				ST1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37		E
				1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2		
				X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DLT		
	Help															
	Cancel (<<< Back														NEXT	
		_												@ 200	0 Terran	

**Difference Data** (Differensdata) väljs automatiskt i rutan **Input data** (Indata). Om du tidigare har bekräftat definitionen av en blankreducering namnger programmet den automatiskt **Blank reduction** (Blankreducering) (se transformerad data i kontrollfältet).

I 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.

För ytterligare detaljer och förklaringar om definitionen och tilldelningen av transformationer, se Magellans användarhandbok.



### 6. Exempel på tillämpning

SM1_9	Prov, försöksgrupp nummer 1, prov-ID nummer 9.
2/2	Antal replikat är 2, totala antal replikat är 2.
x-BL1 eller 1	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 välj **Conc., Dil., Ref.-value** (Koncentration/spädning/referensvärde) från alternativet **Method layout** (Metodlayout) för att definiera respektive värden enligt beskrivningen i provsatsen.

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 från försöksgrupp 1. 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:





## Standardkurva

l kontrollfältet, klicka på **Standard curve** (Standardkurva) från alternativet **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.

Data	Analysis type	ntercepts Axis Graph
	Input data:	Blank reduction
	💿 Standards fro	m Layout
	🔿 Standards fro	m ext. file: Select
	🔿 Standards fro	m experimental group:
	🔿 No Standard	curve graph
	Additional co	oncentrations

På fliken Analysis type (Analystyp), välj linjär regression.

Point to point     Dilinear regression     Non-linear regression     Lubic spline	Data scaling: Lin(x)Lin(y)
<ul> <li>Akima</li> <li>Polynomial</li> <li>Eour parameters</li> <li>Eour parameters Marquardt</li> <li>Five parameters</li> <li>LogitLog</li> </ul>	More
□ Include (0,0)	



## 6. Exempel på tillämpning

På fliken **Axis** (Axel), definierar du märkningen och skalningen av axlarna enligt nedan:

a 🛛 Analysis type	/ Intercepts Axis Graph	
X-axis		
Label:		
Color:	Log. Scaling	
<ul> <li>Auto sel</li> </ul>	lect range	
🔘 Range	Min.: Max.:	
🗹 Grid	Color: Style:	
-Y-axis		
Label:	Blank reduction	
Color:	Log. Scaling	
<ul> <li>Auto sel</li> </ul>	lect range	
🚫 Range	Min.: Max.:	
💌 Grid	Color: Style:	

På fliken **Graph** (Diagram) kan du göra inställningar för diagramrubrik, kurvor, teckensnitt samt diagramvisning.

Label:	IgM-ELISA			
Color:				
Curves				
		Label: Grp. 1		
Color: Symbi Size:	bl: Medium V	Hide curve		
Font		Display		
⊙ Sma		🗹 Legend	Intercepts	
O Med	e	☑ Base points	Error bars	



## Define Cutoffs (Definiera cutoffs)

Välj **Cutoff definition** (Definition av cutoff) i kontrollfältet från alternativet **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/ml	Positiv

Använd följande procedur för att definiera lämpliga gränsvärden:

I rutan Input data (Indata) välj Mean conc. (Medelkoncentration) (UA/ml).

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. Lägg till en färg med hjälp av den nedrullningsbara färgpaletten

- Positiv Röd
- Gränsfall– Blå
- Negativ Grön

Skärmen innehåller följande:

sitive	22		
ermedial	22		
	18	 	
egative			
<b></b>			
<b></b> _			
<b></b>			
<b></b>			
━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━━			
	gative		

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 alternativet **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 första raden skriver du in NC1\_1<8



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:

Validation Con	ditions				
NC1_1<8					
NC1_2<8					
1					
Formula input					
Variable	Operators	Functions			
	00010(010	T GILCUOITS			
BL1		and	~		
BL1 Plate to Plate QC	▼ +	and			
BL1 Plate to Plate QC Input data:	Difference data	and	~		
BL1 Plate to Plate QC Input data:	Difference data	and Mean:	×	1 [	
BL1 Plate to Plate QC Input data: Control 1: Control 2:	Difference data	Mean:			· ]
BL1 Plate to Plate QC	Difference data	Mean:		] <u>†</u>	· _
BL1 Plate to Plate QC Input data: Control 1: Control 2: Control 3:	Difference data	Mean: Mean:	•           •           •           •           •           •           •           •           •           •           •           •           •           •	1 <b>1</b>	·
BL1 Plate to Plate QC Input data: Control 1: Control 2: Control 3: Control 4:	Difference data	Mean: Mean:			
BL1 Plate to Plate QC Input data: Control 1: Control 2: Control 3: Control 4:	Difference data	Mean: Mean: Mean: Mean:			
BL1 Plate to Plate QC Input data: Control 1: Control 2: Control 3: Control 4:	Difference data	Mean: Mean: Mean: Mean:	S:           S:           S:           S:           S:           S:	 	
BL1 Plate to Plate QC Input data: Control 1: Control 2: Control 3: Control 4:	Difference data	Mean:			



# 6.2.3 Organisera rapport

Klicka på **Printed report** (Utskriven rapport) i kontrollfältet från alternativet **Data handling** (Datahantering). Den följande skärmen visas:

List      Matrix	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.
vailable data:	Selected data:
	Insert Append Ap

På fliken **Data selection** (Dataurval), finns alla data som är tillgängliga för rapporten i rutan **Available data** (Tillgänglig data). Med hjälp av knapparna **Insert and Append** (Lägg till 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 datan 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 du skapar rapporten måste det förinställda Vertical list/Difference data (Vertikal lista/differensdata) tas bort från rutan Selected data (Utvalda data). Därefter återstår endast Measurement parameters (Mätparameterar) i rutan Selected data (Vald data). Print as List (Skriv ut som lista) måste ändras till Print as Matrix (Skriv ut som matris).



#### 6. Exempel på tillämpning

Print as List   Matrix	Pick an item fror Droj	You can use drag & drop to create a report. n the available data list and drag it into the s o it into the appropriate position inside the rep	elected data list. port.
vailable data:	Si	elected data:	
Instrument data     Reduced data     Difference data     Difference data     Difference data - Mean     Difference data - Standard de     Difference data - Variation co     Transformed data     Concentrations     Qualitative Results     Audit Pasult     Measurement parameters     Remarks     Error protocol     E Audit Trail	Insert Append	E Data Collection	Up Down Properties Remove all
	Import		

Välj **Method layout** (Metodlayout)/**Layout** i rutan **Available data** (Tillgänglig data) och bifoga den som en matris till rapporten genom att klicka på **Append** (Bifoga). Lägg därefter till **Blank reduction** (Blankreducering), **Mean conc.** (Medelkoncentration) **(UA/mI)** samt **Cutoff results** (Cutoff-resultat) i denna matris genom att välja de motsvarande alternativen och sedan klicka på **Insert** (Lägg till).

Append (Bifoga) Graph (Diagram): Standard curve (Standardkurva), Cutoff definition (Definition av cutoff), och QC Validation criteria (QC-kriterier för validering) till den utvalda datan. 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) kan du definiera layouten på rapportens sidhuvud och sidfot (se Magellans användarhandbok för ytterligare detaljer).





## Data Export (Dataexport)

Klicka på **Data export** (Dataexport) i kontrollfältet från alternativet **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** (Tillgänglig data). Klicka på  $\rightarrow$  pilen för att lägga till dem i fönstret **Selected data** (Utvalda data). Skärmen visar följande information:

Available data:			Selected data:	
🕀 🛍 Instrument data	>		Layout	Up
🖨 🤆 Reduced data				Down
📃 Difference data				Domi
Difference data - Mean				
Difference data - Standard deviation				
Difference data - Variation coefficie				
🕀 🕰 Transformed data				
🖶 🗍 Concentrations		•		
Cutoff results				
🗊 🖓 'z, Sample IDs		(·		
🖮 🛲 Method layout				
Well positions				
Strip method names				
Original Concentrations				
Dilution factors	v			
Export Options Export to ASCII File		Export to	o Excel	



Anmärkning Data som exporteras bör alltid innehålla layout eller prov-ID.

## Automatiserad datahantering

Klicka på **Automated Data Handling** (Automatiserad datahantering) i kontrollfältet från alternativet **Data handling** (Datahantering).

Mara
More
More
More

Välj export to ASCII file (exportera till ASCII-fil), och view results after measurements (visa resultat efter mätningar). I Magellan Tracker, väljs save workspace (spara arbetsyta) 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/Ed	lit a Method				$\mathbf{X}$	
Save in:	The meth	Name 🔺	Remarks	Status		
Filename:	Method1.mth					
File remark:	s:			~		
				<u>~</u>		
Audit trail c	omment:					
Org	anize favorites Audit trail	Signatures	Method password:			
H Ca	telp ancel <<< Back	🗌 Run this	method now SAVE&	FINISH		
			© 20	109 Tecan		
Filena textfält	t <b>me</b> (Filnamn)	Ett filnamn måste a föreslås automatisk	nges. Ett stand t, men detta ka	dardfilnamn an ändras.		
File re (Filkon textfält	e <b>marks</b> nmentarer) t	De kommentarer so och visas tillsamma	om anges här k Ins med filnam	kommer att spa net.	aras	
Audit komm textfält	trail nentar t	De kommentarer so i Audit trail. Detta a <b>Magellan Tracker</b> .	om anges här k Iternativ är bar	kommer att spa a tillgängligt i	aras	
<b>Organize Favorites</b> (Organisera favoriter) knapp		Dialogrutan <b>Organize Favorites</b> (Organisera favoriter) visas.				
Metho (Löser textfält	od password nord för metod) t	Ange ett lösenord fo	ör att skydda n	netoden.		
Run th now (l	h <b>is method</b> Kör denna metoden vssruta	Metoden kommer a på Save&Finish (Sp	tt köras efter a bara och slutfö	tt du har klicka r).	at	
nu) kij						

För ytterligare detaljer, se Magellans användarhandbok.



## 6.2.4 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). I dialogen **Start Measurement** (Starta mätning) tilldelas ett standardnamn på arbetsytan och kan ändras om så önskas.

Start Measurement				×
Measurement			Measurement parameters	
Workspace: Method: Sample ID list: Instrument Use sta Plate in Movements Please note:	05112009-001.wsp Method1.mth cker Temp. control Injector control	Arb. cycle kin Modify layout Insert Current: n. def °C Target: n. def °C Optimize Z-position	Plate Plate Description: [GRE96it] - Greiner 96 Flat Tr Plate with Cover: No Barcode: No Patt of Plate Range: A1:H12 Absorbance Measurement wavelength: 492 nm Reference wavelength: 620 nm Label: Label1	
Help Cancel	<<< Back		START	
			© 2009 Tecan	

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. Felmeddelanden kan förekomma när du utför en mätning utan motsvarande vätskor (t.ex. standarder).



# 6.2.5 Utvärdera resultat

Välj **Evaluate results** (Utvärdera resultat) för att se och utvärdera rådata. Du kan se utvärderingsparametrarna och utvärdera data på nytt.

Detta avsnitt leder dig igenom guiden **Evaluate Results** (Utvärdera resultat) med hjälp av ett exempel på arbetsyta som installerats automatiskt vid installation av Magellan.

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\_InfiniteF50.wsp** från fillistan och klicka på **Make your selection** (Gör ditt val). Beräkningarna utförs och följande plattlayout visas:

Evaluate results - Quantitative BLISA example_Sourise_InfiniteF50.wsp - Quantitative BLISA example_Sourise.mth File Edit Indoument															
e edt		10% 100	1% Zoor	n											
(															
Measurement data 🛞 Measurement data Reference data	Evaluate n		1	2	3	4	5	6	7	8	9	10	11	12	
Measurement data - Mean s - Measurement data v - Measurement data Reference data - Mean	esults	A	BL1 1/1 0.004	ST1_3 2/2 0.207	SM1_1 2/2 0.1	SM1_5 2/2 0.816	SM1_9 2/2 0.174	SM1_13 2/2 0.166	SM1_17 2/2 0.083	SM1_21 2/2 0.085	SM1_25 2/2 0.085	SM1_29 2/2 0.08	SM1_33 2/2 0.162	SM1_37 2/2 0.131	
s - Reference data v - Reference data Measurement data as Colors Reference data as Colors	Edit method	B	NC1 1/2 0.069	ST1_4 1/2 0.418	SM1_2 1/2 0.784	SM1_6 1/2 0.212	SM1_10 1/2 0.196	SM1_14 1/2 0.156	SM1_18 1/2 0.106	SM1_22 1/2 0.216	SM1_26 1/2 0.123	SM1_30 1/2 0.103	SM1_34 1/2 0.131	SM1_38 1/2 0.152	
Reduced data		C	NC1 2/2 0.068	ST1_4 2/2 0.418	SM1_2 2/2 0.764	SM1_6 2/2 0.205	SM1_10 2/2 0.193	SM1_14 2/2 0.155	SM1_18 2/2 0.104	SM1_22 2/2 0.21	SM1_26 2/2 0.129	SM1_30 2/2 0.112	SM1_34 2/2 0.125	SM1_38 2/2 0.155	
Difference data - Mean s - Difference data v - Difference data Difference data as Colors		D	ST1_1 1/2 0.052	ST1_5 1/2 0.838	SM1_3 1/2 0.64	SM1_7 1/2 0.083	SM1_11 1/2 0.098	SM1_15 1/2 0.108	SM1_19 1/2 0.121	SM1_23 1/2 0.11	SM1_27 1/2 0.162	SM1_31 1/2 0.105	SM1_35 1/2 0.127	SM1_39 1/2 0.093	
Transformed data 😒		Ε	ST1_1 2/2 0.051	ST1_5 2/2 0.84	SM1_3 2/2 0.629	SM1_7 2/2 0.085	SM1_11 2/2 0.1	SM1_15 2/2 0.11	SM1_19 2/2 0.125	SM1_23 2/2 0.115	SM1_27 2/2 0.166	SM1_31 2/2 0.111	SM1_35 2/2 0.129	SM1_39 2/2 0.09	
Qualitative Results 😒		F	ST1_2 1/2 0.103	ST1_6 1/2 1.658	SM1_4 1/2 0.323	SM1_8 1/2 0.104	SM1_12 1/2 0.078	SM1_16 1/2 0.153	SM1_20 1/2 0.143	SM1_24 1/2 0.165	SM1_28 1/2 0.112	SM1_32 1/2 0.094	SM1_36 1/2 0.135	SM1_40 1/2 0.143	
QC Validation 🛞 QC Validation criteria		G	ST1_2 2/2 0.103	ST1_6 2/2 1.655	SM1_4 2/2 0.314	SM1_8 2/2 0.099	SM1_12 2/2 0.079	SM1_16 2/2 0.155	SM1_20 2/2 0.136	SM1_24 2/2 0.164	SM1_28 2/2 0.116	SM1_32 2/2 0.092	SM1_36 2/2 0.124	SM1_40 2/2 0.149	AN Austria
Miscellaneous 🛞 Remarks Error protocol Audit Trail Signature		H	ST1_3 1/2 0.206	SM1_1 1/2 0.105	SM1_5 1/2 0.845	SM1_9 1/2 0.199	SM1_13 1/2 0.167	SM1_17 1/2 0.085	SM1_21 1/2 0.081	SM1_25 1/2 0.082	SM1_29 1/2 0.081	SM1_33 1/2 0.167	SM1_37 1/2 0.127		TEC
Meaparant parameters and a second sec															
Heb Carcel (CBack NEXT)															

I varje enskild brunn visas det beräknade värdet. Plattlayouten ändras på motsvarande sätt, beroende på det valda alternativet i kontrollfältet. Parametrar och inställningar kan ändras med hjälp av alternativen 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	_10 SM1_14 SM1_18					
2 0.1	Summary Details Edit					
SM1 1 0.0	Edit kinetic settings Copy kinetic settings Paste kinetic settings					
SM <sup>2</sup> 2 0	Graph: Multilabel Graph: Kinetics Graph: Multilabel kinetics Graph: FLT curves					
SM <sup>r</sup>	Graph: Spectra Graph: Dilution series					
0.0	Mask/Unmask selection Show/Hide layout					

Om du väljer **Summary** (Sammanfattning) visas följande fönster med detaljerad information om definitionen och inställningarna för den valda brunnen:

Well: C5	$\mathbf{X}$						
📮 👜 Method layout							
🕀 🛲 Identifier							
🕀 👜 Liquid Alias							
🕀 🛲 Replicate							
🗉 🛲 Multilabel data reduction	🛨 📠 Multilabel data reduction						
🛨 📖 Transformation formula(s)							
Kinetic transformation formula(s)							
Concentration transformation formula(s)							
🕀 📠 Strip method names							
🗉 🛄 Instrument data							
🗉 🔆 Reduced data							
🗄 🔟 Averages							
🕀 🔏 Transformed data							
🗄 🗊 Concentrations							
💼 🛕 Statistics							
left Up right							
Expand All	ОК						

Klicka på **Next** (Nästa) i fönstret för plattlayouten och dialogrutan **Save as** (Spara som) visas, där kan du 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 **Finish** (Slutför) till höger längst ner på skärmen för att spara filen och stänga guiden. Programmet går tillbaka till guide-listan.



# 6.2.6 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) (Koncentration/spädning/referensvärde)

Utvald identifierare: ST

Enhet: UA/ml

ST1_1	5	(ST1_1Standard 1 första försöksgrupp)
ST1_2	10	(ST1_2Standard 2 första försöksgrupp)
ST1_3	20	(ST1_3Standard 3 första försöksgrupp)
ST1_4	40	(ST1_4Standard 4 första försöksgrupp)
ST1_5	80	(ST1_5Standard 5 första försöksgrupp)
ST1_6	160	(ST1_6Standard 6 första försöksgrupp)

#### 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: Medelkoncentration. (UA/ml)

Begränsningar: 22

18

Positiv ≥ 22 > gränsfall ≥ 18 > negativ

Icke-kompetitivt test



# 5. QC Validation (QC-validering)

Definitioner i Magellan (Kontrollfält – Evaluate data/QC validation) (Utvärdera data/QC-validering):

Indata: Enstaka koncentration (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



# 7. Rengöring, underhåll, och avfall

# 7.1 Inledning

Detta kapitel innehåller följande procedurer:

- Rengöring av instrumentet
- Desinficering av instrumentet
- Underhållning av instrumentet
- Insättning eller ersättning av filter i filterhjulet
- Avfallshantering



# 7.2 Rengöring av instrumentet



#### VARNING RENGÖRINGSPROCEDUREN SKA UTFÖRAS I EN VÄLVENTILERAD LOKAL AV BEHÖRIG OCH UTBILDAD PERSONAL SOM BÄR ENGÅNGSHANDSKAR SAMT

SKYDDSGLASÖGON OCH SKYDDSKLÄDER.

enheten är mycket smutsig, rengör med en trasa fuktad med max 70 % etanol eller ett milt rengöringsmedel, Microcide SQ eller Decon 90. Torka av med en pappersduk.

Om vätska spills på instrumentet ska den omedelbart tas bort för att förhindra att vätska rinner in i det optiska systemet och orsakar prestandaförlust eller ett felmeddelande.



# 7.3 Desinfektion av instrument



#### VARNING

#### OM VÄTSKA SOM SPILLTS PÅ PLATTHÅLLAREN ÄR POTENTIELLT SMITTFÖRANDE SKA DEN DESINFICERAS I ENLIGHET MED GÄLLANDE NATIONELLA LAGAR OCH REGELVERK.

Alla delar av instrumentet som har kommit i kontakt med biologiska prover, patientprover, positiva kontrollprover eller farligt material måste behandlas som potentiellt smittförande.



#### VARNING

DESINFEKTIONSPROCEDUREN OCH DESINFEKTIONSMEDLEN SKA ÖVERENSSTÄMMA MED GÄLLANDE NATIONELLA LAGAR OCH REGELVERK.



#### VARNING

DET ÄR MYCKET VIKTIGT ATT INSTRUMENTET DESINFICERAS NOGGRANT INNAN DET LÄMNAR LABORATORIET ELLER INNAN NÅGON SERVICE UTFÖRS PÅ DET.

Innan instrumentet returneras till den lokala försäljningsrepresentanten eller till ett servicecenter måste alla ytor och platthållaren desinficeras, och ett säkerhetsintyg måste fyllas i av den verksamhetsansvariga. Om ett säkerhetsintyg inte tillhandahålls kan det hända att instrumentet inte godkänns av den lokala säljaren eller servicecentret, eller att tullmyndigheterna behåller det.

# 7.3.1 Desinfektionslösningar

Instrumentets ytor samt instrumentets platthållare ska desinficeras med hjälp av en desinficeringslösning som:

- Microcide SQ
- Decon 90
- 70 % etanol



# RISK FÖR BRAND OCH EXPLOSION!

VARNING

ALKHOL, SOM ETANOL ELLER ISOPROPANOL ÄR BRANDFARLIGT OCH KAN ORSAKA EXPLOSIONER OCH/ELLER BRAND VID FELAKTIG HANTERING. VIDTA LÄMPLIGA FÖRSIKTIGHETSÅTGÄRDER FÖR LABORATORIER.



FÖRSIKTIGHET ANVÄND ALDRIG ACETION EFTERSOM DET SKADAR HÖLJENA.



# 7.3.2 Desinfektionsprocedur

Om laboratoriet saknar specifika desinfektionsprocedur ska följande procedur användas för desinfektion av utsidan och platthållaren på instrumentet.







### VARNING DESINFEKTIONSPROCEDUREN SKA UTFÖRAS I EN VÄLVENTILERAD LOKAL AV BEHÖRIG OCH UTBILDAD PERSONAL SOM BÄR ENGÅNGSHANDSKAR SAMT SKYDDSGLASÖGON OCH SKYDDSKLÄDER.

#### FÖRSIKTIGHET YTDESINFICERINGSMEDLET KAN PÅVERKA DITT INSTRUMENTS PRESTANDA NEGATIVT OM DET TILLÄMPAS PÅ INSTRUMENTETS INSIDA ELLER OAVSIKTLIGT TRÄNGER IN I INSTRUMENT.

#### VARNING

KOPPLA BORT INSTRUMENT FRÅN STRÖMFÖRSÖRJNINGEN FÖRE DESINFICERING FÖR ATT UNDVIKA RISK FÖR BRAND ELLER EXPLOSION.

- 1. Använd skyddshandskar, skyddsglasögon och skyddskläder.
- 2. Förbered en lämplig behållare för alla engångsartiklar som används under desinfektionsproceduren.
- 3. Koppla ur instrumentet från programvaran och stäng av instrumentet.
- 4. Flytta försiktigt platthållaren ut ur instrumentet.



- 5. Applicera försiktigt desinficeringslösning enligt tillverkarens användarhandbok på platthållaren. Använd inte för mycket desinfektionsmedel för att förhindra att lösningen flödar in i instrumentet eller smutsar ner linserna när du flyttar in platthållaren in i apparaten.
- 6. Efter nödvändig kontakttid (enligt tillverkarens användarhandbok) torka av platthållaren med en mjuk pappershandduk fuktad med ett milt rengöringsmedel eller destillerat vatten för att ta bort alla spår av desinfektionsmedlet.



7. Placera försiktigt platthållaren in i instrumentet.



- 8. Applicera försiktigt desinfektionsmedel enligt tillverkarens användarhandbok på platthållarens basplatta.
- 9. Efter nödvändig kontakttid torka av platthållarens basplatta med en mjuk pappershandduk fuktad med ett milt rengöringsmedel eller destillerat vatten för att ta bort alla spår av desinfektionsmedlet.
- 10.Applicera försiktigt desinficeringslösning enligt tillverkarens användarhandbok på alla ytor på utsidan av instrumentet.
- 11.Efter nödvändig kontakttid torka av instrumentet med en mjuk pappershandduk fuktad med ett milt rengöringsmedel eller destillerat vatten för att ta bort alla spår av desinfektionsmedlet.
- 12.Torka av ytorna på utsidan av instrumenten torra med en mjuk pappershandduk.
- 13.Upprepa desinfektionsproceduren på alla tillbehör som flyttas eller skickas tillbaka.
- 14.Kasta behållaren med engångsartiklar i enlighet med gällande nationella lagar och regelverk.
- 15.Desinficera dina händer och tvätta dem med ett milt tvättmedel.

När du skickar tillbaka instrumentet till den lokala säljaren/servicecentret fortsätter du med följande steg:

16.Packa in instrumentet och dess tillbehör.

17.Fyll i säkerhetsintyg (se nedan) och fäst den på utsidan av lådan så att den är tydligt synlig.

## 7.3.3 Säkerhetsintyg

För att garantera personalens säkerhet och hälsa ombeds våra kunder att fylla i ett **Säkerhetsintyg** (som levererades tillsammans med instrumentet) och fästa en kopia på behållaren där instrumentet returneras (synligt från utsidan av transportbehållaren!) och en annan kopia i fraktdokumenten innan instrumentet skickas till servicecentret för service eller reparation.

Instrumentet måste desinficeras på plats innan det fraktas i väg (se 7.3.2 Desinfektionsprocedur).

Desinfektionen måste utföras i ett välventilerat rum av behörig och utbildad personal som använder puderfria engångshandskar, skyddsglasögon och skyddskläder.



Desinfektionen ska utföras i enlighet med nationella, regionala och lokala regelverk.

Om ett säkerhetsintyg inte tillhandahålls kan det hända att instrumentet inte godkänns av servicecentret.

Din lokala Tecan kundsupport kan vid behov skicka en ny kopia av säkerhetsintyget till dig.

# 7.4 Förebyggande underhållschema för INFINITE F50 PLUS

Följande förebyggande underhållsprocedur rekommenderas.

# 7.4.1 Månatlig

Rengör höljet och platthållaren med ett milt rengöringsmedel åtminstone en gång per månad, oftare vid behov.



## FÖRSIKTIGHET ANVÄND ALDRIG ACETON EFTERSOM DET SKADAR HÖLJENA.

# 7.4.2 Vart fjärde år

Det rekommenderas att filtren byts ut vart fjärde år.

# 7.5 Filterbyte och installation

INFINITE F50 PLUS måste vara ansluten till Magellan för att kunna göra en filterinsättning/filterbyte guidad av en programvara. Om anslutningen förloras under proceduren på grund av att anslutningen mellan instrumentet och datorn avbryts oavsiktligt måste Magellan avslutas och instrumentet måste stängas av. Fortsätt i så fall med proceduren enligt nedan. Återupprätta du anslutningen när du är klar genom att starta om enheten och Magellan, och definierar de nyinsatta filtren.



## FÖRSIKTIGHET

VAR FÖRSIKTIG NÄR DU HANTERAR FILTRENA SÅ ATT DE INTE BLIR REPADE ELLER SMUTSIGA AV FINGERAVTRYCK ELLER DAMM.



## 7.5.1 Steg för filterutbyte

Filtren i standardfilterhjulet kan bytas ut eller kompletteras enligt följande procedure:

- 1. I fönstret Wizard list (Guide-lista), klicka på Miscellaneous (Övrigt).
- 2. Klicka på Instrument control (Instrumentetkontroll)
- 3. Klicka på Define filter slides... (Definiera filterhållare)
- 4. Klicka på Filter switching (Filterutbyte) för att starta proceduren.
- 5. Ta bort eventuella mikroplattor från platthållaren!
- Luta instrumentet försiktigt bakåt tills det ligger på baksidan med botten riktad mot dig.
- 7. Ta bort plattskyddet från instrumentets botten genom att ta bort de orangea skruvarna.



8. Ta bort det magnetiskt fastsatta filterhjulet genom att försiktigt dra ut det ur instrumentet.



9. Placera filterhjulet på en ren plan yta.



10.När du byter ut ett filter, använd verktyget för filtersammansättning för att ta bort filtret från filterspåret.

Kontakta din lokala säljare för Tecan för verktyget för filtersammansättning och tillgängliga filter.



- 11. Rikta verktyget för filtersammansättning med stoppringens skåra. Vrid verktyget och ta bort stoppringen genom att dra ut den ur filterspåret.
- 12.Vänd filterhjulet så att filtret glider ut ur spåret. Använd inte verktyget för filtersammansättningen för att trycka ut filter ur filterspåret, eftersom filtret kan bli förstöras.
- 13.Ett nytt filter måste sättas in i filterspåret i korrekt riktning. Se till att inte skrapa filtret eller lämna fingeravtryck på det.





Anmärkning Kontrollera att filtret är korrekt isatt.



14.Placera stoppringen på ändan av verktyget för filtersammansättning och vrid på den så att den inte kan glida av.



- 15. Använd verktyget för filtersammansättning för att trycka in stoppringen i filterspåret och tryck den stadigt på plats.
- 16.Rotera verktyget tills skåran i stoppringen är i linje med ändan på verktyget för filtersammansättning och ta bort verktyget.



- 17.Placera filterhjulet tillbaka i filterhjulspåret och tryck in det djupare tills det är magnetiskt fastsatt.
- 18.Sätt tillbaka plattskyddet på instrumentets botten med de tre orangea skruvarna.
- 19. Lägg enheten tillbaka till ett uppåtvänt läge.
- 20.Klicka på OK för att avsluta proceduren och initialisera filterhjulet.
- 21. Definiera det nya filtret (se nästa kapitel för detaljerad procedur).


#### 7.5.2 Definiera filter

I dialogrutan **Filter Definition** (Filterdefinition) ska du tilldela de ersatta filtren lämpliga våglängder genom att ange de nya våglängderna i motsvarande positioner.

Om ett filter har satts in i en ny filterposition aktiverar du lämplig filterposition genom att markera kryssrutan och ange lämplig våglängd.

Genom att klicka på **Save** (Spara) sparas filterdefinitionerna, och filtren initialiseras.

När filtren har initialiserats är instrumentet redo för mätningar.



Anmärkning Se till att inte förväxla filterpositioner och filtervåglängder, eftersom detta leder till felaktiga mätdata.

# 7.6 Avfall

### 7.6.1 Inledning

Följ laboratorierutiner för hantering av biologiskt avfall enligt nationella eller lokala regelverk.

Detta kapitel tillhandahåller instruktioner om korrekt hantering av förbrukat material som uppstår vid användningen av INFINITE F50 PLUS.



#### FÖRSIKTIGHET FÖLJ ALLA STATLIGA, REGIONALA OCH LOKALA MILJÖBESTÄMMELSER.

#### 7.6.2 Avfallshantering av förpackningsmaterial

Förpackningsmaterialet består av återvinningsbart material. Om du inte har för avsikt att behålla det för framtida användning, till exempel för transport- och förvaringsändamål, kassera förpackningsmaterialet i enlighet med lokala regelverk.

## 7.6.3 Avfallshantering av förbrukat material

#### VARNING

BIOLOGISKA FAROR KAN VARA FÖRKNIPPADE MED AVFALLSMATERIALET (MIKROPLATTOR) SOM UPPSTÅR VID PROCESSER SOM KÖRS PÅ ABSORBANSLÄSAREN INFINITE F50 PLUS.



BEHANDLA DEN ANVÄNDA MIKROPLATTAN SAMT ALLA ENGÅNGSARTIKLAR OCH ALLA SUBSTANSER SOM INGÅTT I PROCESSEN I ENLIGHET MED RIKTLINJERNA FÖR GOD LABORATORIEPRAXIS.

TA REDA PÅ LÄMPLIGA INSAMLINGSPLATSER OCH VILKA METODER FÖR AVFALLSHANTERING SOM ÄR GODKÄNDA LOKALT.



#### 7.6.4 Avfallshantering av instrumentet

Kontakta Tecans lokala kundsupport om du har några frågor om avfallshantering

Föroreningsgrad	2 (IEC/EN 61010-1)
Avfallshantering	Kontaminerat avfall

#### **OBSERVERA**

DIREKTIV 2012/19/EU OM AVFALL SOM UTGÖRS AV ELLER INNEHÅLLER ELEKTRISKA ELLER ELEKTRONISKA PRODUKTER (WEEE)

RISK FÖR ICKE ÖNSKVÄRDA MILJÖEFFEKTER TILL FÖLJD AV AVFALLSHANTERING.

- KASTA INTE ELEKTRISK OCH ELEKTRONISK UTRUSTNING BLAND VANLIGT HUSHÅLLSAVFALL.
- UTTJÄNT ELEKTRISK OCH ELEKTRONISK UTRUSTNING SKA SAMLAS IN SEPARAT.

#### VARNING

BEROENDE PÅ TYP AV TILLÄMPNING KAN DELAR AV INFINITE F50 PLUS HA KOMMIT I KONTAKT MED BIOLOGISKT RISKMATERIAL.

- SE TILL ATT HANTERA DETTA MATERIAL ENLIGT GÄLLANDE SÄKERHETSSTANDARDER OCH BESTÄMMELSER.
- FÖRE AVFALLSHANTERINGEN SKA ALLTID SAMTLIGA DELAR DEKONTAMINERAS.

För mer information om produkten, kontakta:

Tecan Austria GmbH

Untersbergstrasse 1A

A-5082 Grödig/Salzburg

ÖSTERRIKE/EUROPA

T +43 6246 8933 444

E-post: <u>expertline-at@tecan.com</u>

www.tecan.com







# 8. Felsökning

# 8.1 Inledning

Den interna mikroprocessorn styr och kontrollerar elektroniska funktioner samt mätningar, drift och resultat. Om mikroprocessorn upptäcker ett fel eller ett felaktigt tillvägagångssätt visas ett felmeddelande på datorn.

# 8.1.1 Tabell med felmeddelanden och felsökning

I följande tabell ges en kortfattad beskrivning av felmeddelandena och felsökningsåtgärderna.



Anmärkning Om andra felmeddelanden än de som visas i tabellen nedan förekommer skall du kontakta din lokala Tecan kundsupport.

Felmeddelande	Beskrivning	Felsökning
Systemfel		
"Lid Open Error"	Locket är öppet i början av en mätningen	Stäng locket och påbörja mätningen igen
"MTP Init Error"	MTP-transport kunde inte initialiseras	Maskinvaruproblem: Elektronisk defekt, remmen är trasig, MTP- transporten kan vara mekaniskt blockerad.
"MTP lost steps abs( <i>steploss</i> ) > <i>max_steploss</i> "	MTP tappade steg under mätningen Steploss: antal tappade steg max_steploss: antal tillåtna steg	Maskinvaruproblem: Elektronisk defekt, ojämn mekanik
"Filter lost steps abs( <i>steploss</i> ) > <i>max_steploss</i> "	Filterhjulet tappade steg under mätningen. Steploss: antal tappade steg max_steploss: antal tillåtna steg	Maskinvaruproblem: Elektronisk defekt, ojämn mekanik
"USB timeout"	Timeout för USB-kommunikation	Systemfel - rapportera till kundsupport
"Lamp Low! Minimum: <i>minimum</i> , Maximum: <i>maximum</i> " Detta meddelande visas med versioner av fasta programvarurversioner upp till V1.11.	Den uppmätta ljusintensiteten nådde inte upp till det förväntade intervallet mellan <i>minimum</i> och <i>maximum</i>	Maskinvaruproblem: Elektronisk defekt, trasig fiber
I version V1.12 av den fasta programvaruversionen och högre ersätts meddelandet "Lamp Low" med: "Prepare REF check"	Den uppmätta ljusintensiteten nådde inte upp till det förväntade intervallet mellan <i>minimum</i> och <i>maximum</i>	Maskinvaruproblem: Elektronisk defekt, trasig fiber
"Wavelength Not Available! Wavelength: <i>wavelength</i> nm"	Filtret med våglängden <i>wavelength</i> kunde inte hittas i filterhjulet.	Systemfel - rapportera till kundsupport



#### 8. Felsökning

Felmeddelande	Beskrivning	Felsökning
Systemfel		
"Channel Low! Channel: <i>channel_nr</i> , Minimum: <i>minimum</i> , Maximum <i>maximum</i>	Signalen på kanal <i>channel_nr</i> nådde inte det förväntade området mellan <i>minimum</i> och <i>maximum</i>	Maskinvaruproblem: Elektronisk defekt, trasig fiber
"Invalid Wavelength! Wavelength: wavelengthnm"	Filtrets våglängd ligger utanför våglängdsområdet för vit och blå LED	Systemfel - rapportera till kundsupport
"Lamp Overflow!" Minimum: <i>minimum</i> , Maximum: <i>maximum</i> "	Signalen på ADC överskrider det förväntade området mellan <i>minimum</i> och <i>maximum</i>	Maskinvaruproblem: Elektronisk defekt
"Value Not Set: <i>value -1</i> "	Värdet value är inte inställt.	Systemfel - rapportera till kundsupport
"Filter Init Error"	Filtrertransport kunde inte initialiseras	Maskinvaruproblem: Elektronisk defekt, filterhjulstransport mekaniskt blockerad.

# 8.1.2 Definition av 'Overflow (Överflöde)'

Om resultatet av absorbansmätningen inte ligger inom instrumentspecifikationerna (> 4,0 OD) uppstår ett överflöde och det uppmätta OD-värdet för respektive brunn ersätts med "Overflow" (Överflöde). Detta görs av programvaran och inte av själva instrumentet.

#### 8.1.3 Strömavbrott

Vid strömavbrott händer följande:

Strömavbrott i instrumentet, men inte i styrdatorn (t.ex. dator ansluten till ett system för avbrottsfri strömförsörjning): USB-anslutningen mellan instrumentet och datorn kommer att förloras. Felmeddelande av kontrollprogramvaran skapas.

Strömavbrott sker för både instrumentet och styrdatorn: starta om datorn. Inga mätdata kommer att finnas tillgängliga.

# 9. Förkortningar, varumärken och symboler

# 9.1 Förkortningar

Följande förkortningar tillhandahålls som referens och kan förekomma i användarhandboken.

A	Ampere
AC	Växelström
ADC	Analog-till-digitalomvandlare
ANSI/SBS	American National Standards Institute/Society for Biomolecular Screening
ASCII	American Standard Code for Information Interchange
ASTM	American Society for Testing and Material
°C	grader Celsius
CE	CE-märkning om överensstämmelse
CFR	Code of Federal Regulations (Federala förordningar)
cm	centimeter
DC	Likström
EC	Europeiska gemenskapen
ELISA	Enzymkopplad immunadsorberande analys (Enzyme-Linked Immuno Sorbent Assay)
EN	Europeisk norm
°F	grader Fahrenheit
FDA	Food and Drug Administration (Amerikanska läkemedelsverket)
Hz	Hertz
IEC	Internationella elektrotekniska kommissionen
ID	Identifikation
IFU	Användarhandbok
IQ	Installationskvalificering
IVD	In vitro-diagnostik
IVDR	Förordning (EU) 2017/746 om medicintekniska produkter för <i>in vitro</i> -diagnostik (IVDR)
kg	kilogram
I	liter
LED	Lysdiod
LIS	Laboratorieinformationssystem
mg	milligram
ml	milliliter
mm	millimeter
MTP	mikroplatta
μΙ	mikroliter
NFM	Icke-järnmetaller
NIST	National Institute of Standards and Technology



#### 9. Förkortningar, varumärken och symboler

nanometer
Nationally Recognized Testing Laboratory
Optisk densitet
Driftskvalificering
Tryckt kretskort
Radiofrekvens
Restriction of the Use of Certain Hazardous Substances (Minskning av användningen av vissa farliga ämnen)
Standard Operating Procedure (Standardiserade tillvägagångssätt)
Universal Serial Bus
Godtyckliga enheter
Technischer Überwachungsverein (Provningsanstalt)
Volt
Voltampere
Waste electrical and electronic equipment (Uttjänt elektrisk och elektronisk utrustning)

# 9.2 Varumärken

Följande produktnamn och alla registrerade och oregistrerade varumärken som nämns i det här dokumentet används endast i identifieringssyfte och är respektive ägares exklusiva egendom:

- Magellan<sup>™</sup>, Infinite®, MultiCheck<sup>™</sup>, Tecan® och Tecan-logotypen är varumärken som tillhör Tecan Group Ltd., Männedorf, Schweiz
- Windows® and Excel® är registrerade varumärken som tillhör Microsoft Corporation, Redmond, WA, USA
- Pentium<sup>®</sup> och Atom<sup>™</sup> är registrerade varumärken som tillhör Intel Corporation, Santa Clara, CA, USA
- Adobe® Reader® är ett registrerat varumärke som tillhör Adobe Systems Incorporated, Seattle, WA, USA
- Microcide SQ<sup>TM</sup> är ett registrerat varumärke som tillhör Global Biotechnologies Inc., Portland, ME, USA
- Decon 90<sup>™</sup> är ett registrerat varumärke som tillhör Decon Laboratories Ltd., Hove, East Sussex, Storbritannien



# 9.3 Symboler

Följande symboler visas på instrumentet.

	Tillverkare
$\sim$	Tillverkningsdatum
CE	CE-märkning om överensstämmelse
UK CA	Märkningen United Kingdom Conformity Assessed visar att produkten följer gällande bestämmelser i Storbritannien.
ĺĺ	Se användarhandboken
IVD	Medicinteknisk produkt för in vitro-diagnostik
UDI	Unik enhetsidentifiering UDI-symbolen identifierar databäraren på etiketten.
REF	Katalognummer
SN	Serienummer
●	USB-etikett
X	WEEE-symbol
	Kina RoHS-symbol
c SUD US	NRTL TÜV SÜD MARK
	Biologiska risker



# Register

## Α

Absorbans	23
Användarförfrågan	25
Användarprofil	11
Åtgärder Kontrollfält	23
Avfall	65
Avfallshantering	73
Förbrukat material	73
Förpackningsmaterial	73
Instrument	74

## D

Definiera filter	73
Definiera mätningar	
Del av platta	
Desinfektion	
Lösningar	
Procedur	67
Säkerhetsintyg	
Driftskvalificering	
Linjäritetstest	
Precisionstest	
Tester med mikroplattor	
Tester med MultiCheck	

# Е

Exempel på tillämpning	43
Kvantitativ Elisa	43

## F

75
75
69
35
70
77

## G

Guide	
Bifoga signatur	20
Guide-lista	
Skapa/redigera en metod	
Skapa/redigera en prov-ID-lista	20
Starta mätning	19
Utvärdera resultat	19
Guide-lista	19

#### I

Incubation (Inkubation)	25
Infopanel	26

#### Instrument

Funktioner	
Plats	
Tillbehör	
Instrumentbeskrivning	33

#### Κ

Kinetik Kontrollfält	
Kinetikcykel	
Kinetikförhållande	
Kinetiska mätningar	29
Kommentar	
Kontrollfält	
Åtgärder	
Kinetik	
Laboratorieartiklar	
Mätningar	
Övrigt	
Kontrollfältet	22
Kvalitetskontroll	

#### L

Laboratorieartiklar	Kontrollfält	22

#### Μ

Magellan	. 17
Mätningar Kontrollfält	. 23
Mätningar med multipla inmärkningar	. 28
Mikroplattor	. 37
Optimera prestanda	32
Miliökrav	. 14

#### 0

Optimera prestanda	32
Optimering av prestanda	32
Överflöde	76
Övrigt Kontrollfält	25

#### Ρ

Panel för arbetsflöde	
Platta	22
Programelement	
Absorbans	23
Användarförfrågan	
Del av platta	
Dra in och släppa	
Inkubation	
Kinetikcykel	24
Kinetikförhållande	24
Kommentar	
Platta	
Skakning	23
Vänta - timer	25



## Register

Programvara	
Installationskvalificering	
Programvara	17
Installation	17
Magellan	17
Magellan Tracker	17
Programvara	
Användargränssnitt	
Programvara	
Guide-lista	19

## R

Redigeraren för mätparametrar	. 21
Rengöring	. 65

# S

Säkerhet	7
Säkerhetsintyg	68
Självkontroll	
Śkakning	

Slå PÅ instrumentet	16
Slutpunktsmätningar	27
Specifikationer	36
Strömavbrott	
Strömkrav	14
Symboler	
Systemkrav	15

#### U

Underhåll	
Underhållschema	69
Underhåll	65, 69
Uppacking	
Steg för uppackning	13
Uppackning	
Uppackning och inspektion	13

## V

Vänta - timer 25
------------------



# **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.).

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: Model Designation:	Microplate Reader
	INFINITE F50 PLUS
Article Numbers:	30183570

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

#### Machinery Directive RoHS Directive

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

The Supply of Machinery (Safety) Regulations 2008 The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012

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 Infinite F50 PLUS instruments with the article numbers listed above. The model designation varies depending on the specific model with different article number.