

Manual de utilizare pentru Absorbance Reader

# INFINITE F50 PLUS Software MAGELLAN



Nr. articol document:	30186918
	2023-06
Nr. versiune document:	1.2
Versiune firmware:	3.33
Versiune de software <b>Magel</b> l	l <b>an</b> : 7.5







#### AVERTIZARE CITIȚI CU ATENȚIE ȘI RESPECTAȚI MANUALUL DE UTILIZARE ÎNAINTE DE UTILIZAREA INSTRUMENTULUI.

#### Indicație

S-au depus toate eforturile pentru evitarea erorilor din cadrul textelor şi diagramelor; totuşi, Tecan Austria GmbH nu îşi asumă responsabilitatea pentru orice eroare, care ar putea apărea în acest document.

Conform principiilor sale, Tecan Austria GmbH îşi îmbunătăţeşte produsele, atunci când devin disponibile noi tehnici şi componente. Prin urmare, Tecan Austria GmbH îşi rezervă dreptul de a modifica oricând specificaţiile *în urma verificării, validării şi aprobărilor corespunzătoare.* 

Apreciem orice comentarii referitoare la acest document.



#### Producător

Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria T: +43 6246 89330 www.tecan.com E-mail: office.austria@tecan.com

#### Informații copyright

Conținutul acestui document reprezintă proprietatea a Tecan Austria GmbH și nu trebuie copiat, reprodus sau transferat unei alte persoane sau altor persoane, fără acordul scris în prealabil.

Copyright © Tecan Austria GmbH Toate drepturile rezervate. Tipărit în Austria.

#### Declarația pentru certificatul UE

A se vedea ultima pagină pentru Manualul de utilizare.

#### Despre manualul de utilizare

Instrucțiuni originale. În acest document este vorba de **Manual de utilizare** (IFU) pentru INFINITE F50 PLUS absorbance reader, care este destinat măsurării absorbanței luminoase (densitate optică) a probelor în microplăci cu 96 godeuri. Acestea servesc ca referință și ghid pentru utilizator.

Acest document vă oferă instrucțiuni cu privire la:

- Instalarea instrumentului
- Operarea instrumentului
- Curățarea și întreținerea instrumentului

#### Observații despre screenshot-uri

Numărul versiunii indicat pe screenshot-uri este posibil să nu fie identic cu cel al versiunii curent publicate. Screenshot-urile se înlocuiesc, numai dacă s-a modificat conținutul care are legătură cu aplicația.



#### Istoric de modificări – Instrucțiuni de utilizare

Versiunea	Modificări
V 1.2	Noua DoC (declarație de conformitate)

## Avertizări, precauții și indicații

Următoarele tipuri de indicații sunt utilizate în cadrul acestei publicații pentru a evidenția informațiile importante sau pentru a avertiza utilizatorul cu privire la situațiile potențial periculoase:



## Cuprins

1.	Sigu	ranţa		7
	1.1	Siguranța	a instrumentului	7
2.	Gene	eralităti		9
	2.1	Domeniu	de utilizare	9
	2.2	Principiu	I de functionare	10
	2.3	Profilul u	tilizatorului	11
		2.3.1	Utilizator profesionist - Nivel administrator	11
		2.3.2	Utilizatorul final sau utilizatorul de rutină	11
		2.3.3	Tehnicieni de service	11
3.	Activ	vități preli	minare	.13
	3.1	Despach	etarea si verificarea	13
		3.1.1	Verificarea pachetului livrat	13
		3.1.2	Procedura de despachetare	13
	3.2	Cerințe d	e energie	14
	3.3	Cerințe d	e mediu înconjurător	14
	3.4	Cerințe d	le sistem	15
	3.5	PORNIRE	EA instrumentului	16
	3.6	Software		17
		3.6.1	Introducere / Domeniu de utilizare	17
		3.6.2	Procedura de instalare	17
		3.6.3	Calificarea instalării	18
		3.6.4	Inceperea lucrului cu Magellan	19
	3.7	Magellan	- Editor parametrii de măsură	21
		3.7.1	Control Bar (Bară de control)	22
		3.7.2	Workflow Pane (Panou workflow)	26
		3.7.3	Into Pane (Panou Into)	27
	3.8	Magellan	- Definirea masurarilor	27
		3.8.1	Definirea măsurărilor endpoint	،21 مد
		3.0.Z	Definirea masurarilor multilaber	20 20
		3.0.3 3.8.4	Elemente de program Indeptare si Declansare	29 29
	39	Ontimiza	re nentru nerformantă maximă	
	0.0	391	Amplasarea instrumentului	
		3.9.2	Procedura de operare	32
4	Cara	cteristici	instrument	33
	4 1	Caracteri	istici instrument	.00
		4.1.1	Agitare microplacă	33
	4.2	Descriere	e instrument	33
	4.3	Descriere	e centrifugă cu filtre	35
	4.4	Specifica	ıţii instrument	36
		4.4.1	Specificații generale	36
		4.4.2	Specificații de măsură	37
		4.4.3	Microplăci	37
	4.5	Accesori	i instrument	38
5.	Cont	rol al cali	tății	.39
	5.1	Introduce	ere	39
	5.2	Procedur	a de auto-verificare	39
	5.3	Calificare	e operațională (OQ)	39
		5.3.1	Test MultiCheck	39
		5.3.2	Testarea microplăcii	39
		5.3.3	Precizia instrumentului cu probele de lichid	40
		5.3.4	Linearitatea instrumentului cu probe de lichid	40
6.	Exen	nplu de ap	plicație	.43
	6.1	Introduce	ere	43
	6.2	Exemplu	detaliat: Analiză cantitativă ELISA	43



		6.2.1	Descrierea setului de testare	
		6.2.2	Crearea unei metode	45
		6.2.3	Organizarea raportului tipărit	
		6.2.4	Run the Method (Rulare metodă)	61
		6.2.5	Evaluarea rezultatului	
		6.2.6	Rezumatul definiției ELISA cantitative în Magellan	64
7.	Cură	ţarea, în	ntreținerea și eliminarea deșeurilor	65
	7.1	Introdu	Icere	
	7.2	Curăţar	rea instrumentului	
	7.3	Dezinfe	ectarea instrumentului	
		7.3.1	Soluții de dezinfectare	
		7.3.2	Procedura de dezinfectare	
		7.3.3	Certificat de siguranță	
	7.4	Plan de	e întreținere preventivă pentru INFINITE F50 PLUS	69
		7.4.1	Lunar	
		7.4.2	La fiecare 4 ani	69
	7.5	Înlocuir	rea filtrelor și instalarea	69
		7.5.1	Procedura de schimbare a filtrelor	70
		7.5.2	Definirea filtrelor	73
	7.6	Elimina	area deşeurilor	73
		7.6.1	Introducere	73
		7.6.2	Eliminarea ambalajului	73
		7.6.3	Eliminarea materialelor de lucru	73
		7.6.4	Eliminarea instrumentului	74
8.	Rem	ediere d	lefecțiuni	75
	8.1	Introdu	icere	
		8.1.1	Tabel al mesajelor de eroare și al remedierii defecțiunilor	75
		8.1.2	Definiția 'Overflow'	
		8.1.3	Cădere a tensiunii	
9.	Abre	vieri, ma	ărci înregistrate și simboluri	77
	9.1	Abrevie	eri	
	9.2	Mărci îr	nregistrate	
	9.3	Simbol	 uri	
Index				81
Servi	ciul d	e asiste	nță clienți Tecan	83



# 1. Siguranţa

## 1.1 Siguranța instrumentului

- Respectaţi întotdeauna prescripţiile de bază privind siguranţa, atunci când utilizaţi acest produs pentru a reduce riscul rănirii, incendiului sau şocului electric.
- 2. Citiţi şi înţelegeţi toate informaţiile din Manualul de utilizare (IFU). Dacă nu citiţi, nu înţelegeţi sau nu respectaţi instrucţiunile din acest document, pot rezulta deteriorări ale produsului, rănirea personalului de operare sau performanţa scăzută a instrumentului. Tecan nu poartă responsabilitatea daunelor sau vătămărilor corporale ca urmare a deservirii incorecte a dispozitivului.
- Respectaţi toate indicaţiile de AVERTIZARE şi PRECAUŢIE din acest document.
- 4. Deconectați întotdeauna dispozitivul de la sursa principală de alimentare cu energie electrică înainte de curățarea și dezinfectarea acestuia.
- 5. Nu desfaceți niciodată carcasa instrumentului.
- Respectaţi prescripţiile corespunzătoare de laborator privind siguranţa, cum ar fi purtarea îmbrăcămintei de protecţie (de ex. mănuşi, halat de laborator şi ochelari de protecţie) şi aplicarea procedurilor aprobate de laborator privind siguranţa.



#### PRECAUŢIE

ÎN CAZUL ÎN CARE INSTRUCȚIUNILE INDICATE ÎN ACEASTĂ PUBLICAȚIE NU SUNT RESPECTATE ÎN MOD CORESPUNZĂTOR, INSTRUMENTUL POATE FI DETERIORAT SAU PROCEDURILE NU POT FI EXECUTATE ÎN MOD CORECT ȘI PRIN URMARE, SIGURANȚA INSTRUMENTULUI NU POATE FI GARANTATĂ.

Se presupune că utilizatorii instrumentului dețin experiența necesară și sunt familiarizați cu prescripțiile necesare privind siguranța pentru manipularea chimicalelor și substanțelor biologice periculoase.

Conformitatea cu următoarele legi și directive:

- Legea privind protecția industrială națională
- Reglementările de prevenire a accidentelor
- Fişe tehnice de securitate ale producătorilor de reactivi







ÎN FUNCȚIE DE APLICAȚII, UNELE PIESE ALE DISPOZITIVULUI INFINITE F50 PLUS POT INTRA ÎN CONTACT CU MATERIAL BIOLOGIC PERICULOS/MATERIAL INFECȚIOS.

ASIGURAȚI-VĂ CĂ NUMAI PERSONALUL CALIFICAT UTILIZEAZĂ INSTRUMENTUL. ÎN CAZ DE SERVICE SAU ATUNCI CÂND MUTAȚI SAU UTILIZAȚI INSTRUMENTUL, DEZINFECTAȚI ÎNTOTDEAUNA INSTRUMENTUL ÎN CONFORMITATEA CU INSTRUCȚIUNILE DATE ÎN ACEST DOCUMENT.

RESPECTAȚI PRESCRIPȚIILE DE LABORATOR PRIVIND SIGURANȚA, CUM AR FI PURTAREA ÎMBRĂCĂMINTEI DE PROTECȚIE, ATUNCI CÂND SE LUCREAZĂ CU SUBSTANȚE POTENȚIAL INFECȚIOASE.

#### AVERTIZARE

ACEST INSTRUMENT ESTE ÎN CONFORMITATE CU CERINȚELE PRIVIND EMISIILE ȘI IMUNITATEA DESCRISE ÎN IEC 61326-2-6; TOTUȘI, MEDIUL ELECTROMAGNETIC TREBUIE EVALUAT ÎNAINTE DE UTILIZAREA INSTRUMENTULUI.

ESTE REPONSABILITATEA UTILIZATORULUI DE A SE ASIGURA CĂ SE MENŢINE UN MEDIU ELECTROMAGNETIC COMPATIBIL PENTRU INSTRUMENT, ASTFEL ÎNCÂT ACESTA SĂ FUNCȚIONEZE CONFORM DESTINAȚIEI PREVĂZUTE.

NU UTILIZAȚI INSTRUMENTUL ÎN IMEDIATA APROPIERE A SURSELOR DE RADIAȚII ELECTROMAGNETICE PUTERNICE (DE EX. RF NEECRANATE INTENȚIONAT), DEOARECE ACESTEA POT ÎMPIEDICA FUNCȚIONAREA ADECVATĂ A INSTRUMENTULUI ȘI POT CAUZA DE ASEMENEA REZULTATE INCORECTE.



# 2. Generalități

## 2.1 Domeniu de utilizare

INFINITE F50 PLUS este un cititor automat de absorbanță pentru microplăci cu 96 godeuri care include software-ul Magellan pentru utilizarea profesională în laborator în scopul măsurării absorbanței luminoase (densitate optică) a mediilor lichide omogene pentru utilizarea în diagnosticul in vitro.

Instrumentul este destinat utilizării, în principal, în analiza diagnosticului in vitro a probelor din corpul uman furnizate dintr-un test cu anticorpi marcați enzimatic (ELISA) selectat de utilizator.

Informațiile privind diagnosticul specific și tipul de probă sunt definite de testul selectat.

INFINITE F50 PLUS este destinat măsurării și evaluării analizelor calitative, semicantitative și cantitative, conform parametrilor de diagnosticare prevăzuți în acest sens și specificațiilor instrumentului.

Produsul este destinat utilizării în laborator de către personal instruit. Produsul nu este destinat uzului casnic sau de către nespecialiști.



i



Dacă INFINITE F50 PLUS absorbance reader sau software-ul Magellan a fost modificat în orice măsură, garanția nu va mai fi valabilă și conformitatea regulamentară a instrumentului va fi anulată.

#### Indicație

Autoritatea operațională trebuie să folosească numai seturi de testare cu etichete CE pentru aplicații clinice de diagnosticare. Autoritatea operațională trebuie să se asigure că această combinație a unui set special de testare cu un INFINITE F50 PLUS absorbance reader cu etichetă CE și variantele sale au fost validate pentru a întruni cerințele Regulamentului IVD 2017/746 sau reglementările naționale sau locale corespunzătoare.

Dacă INFINITE F50 PLUS absorbance reader este utilizat pentru diferite aplicații, altele decât cele menționate mai sus sau dacă este utilizat cu alt software decât Magellan, nu mai este conform IVD și autoritatea operațională este responsabilă de utilizarea respectivă și de validarea necesară.



Rezultatele obținute folosind INFINITE F50 PLUS depind de utilizarea adecvată a instrumentului și microplăcilor, conform instrucțiunilor date în acest document, precum și de compușii lichizi utilizați (reactivi, chimie). Instrucțiunile pentru utilizare, depozitare și aplicații, care includ probe sau reactivi, trebuie respectate cu strictețe. Prin urmare, rezultatele trebuie interpretate cu atenție.

Indicație Nu desfaceți niciodată carcasa instrumentului, în caz contrar garanția va fi anulată.







## 2.2 Principiul de funcționare

INFINITE F50 PLUS este un cititor de absorbanță cu 8 canale pentru măsurarea densității optice (OD) a lichidelor din microplăcile cu 96 godeuri.

Lumina creată de sursa de lumină LED traversează un filtru optic din centrifuga cu filtre pentru selectarea lungimii de undă necesare pentru măsurarea absorbanței.

După traversarea filtrului, fasciculul luminos este divizat în opt fibre optice pentru focalizarea pe opt godeuri ale microplăcii.

Fasciculele luminoase care traversează godeurile microplăcii sunt direcționate de opt lentile optice și transmise apoi către opt fotodiode în vederea detecției.

Astfel, cititorul INFINITE F50 PLUS măsoară opt godeuri în paralel, ceea ce corespunde unei coloane ale unei microplăci cu 96 godeuri.

Pentru controlul și reglarea intensității luminoase a lămpii LED este utilizată o fibră optică suplimentară (canal de referință) unde lumina ocolește microplaca.

Valorile de transmisie obținute sunt convertite în valori de densitate optică (OD) conform următoarei formule:

## Transmissi on T = $\frac{I}{I_{o}}$

- Io = intensitatea luminii de incidență
- I = intensitatea luminii transmise

OD este logaritmul transmisiei reciproce.

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



## 2.3 Profilul utilizatorului

### 2.3.1 Utilizator profesionist - Nivel administrator

Administratorul este o persoană care dispune de pregătirea tehnică necesară și de competențe și experiență corespunzătoare. Dacă produsul este utilizat în mod corespunzător, această persoană este capabilă să recunoască și să evite pericolele.

Administratorul dispune de competențe vaste și de capacitatea de instruire a utilizatorului final sau a utilizatorului de rutină în privința protocoalelor analizei în legătură cu un produs Tecan în cadrul domeniului de utilizare.

Sunt necesare competențe de folosire a calculatorului și de limba engleză.

## 2.3.2 Utilizatorul final sau utilizatorul de rutină

Utilizatorul final sau utilizatorul de rutină este o persoană care dispune de pregătirea tehnică necesară și de competențe și experiență corespunzătoare. Dacă produsul este utilizat în mod corespunzător, această persoană este capabilă să recunoască și să evite pericolele. Sunt necesare competențe de folosire a calculatorului și de limbă pentru țara de instalare sau de limba engleză.

## 2.3.3 Tehnicieni de service

Tehnicianul de service este o persoană care dispune de pregătirea tehnică necesară și de competențe și experiență corespunzătoare. Dacă produsul necesită lucrări de service sau de întreținere, această persoană este capabilă să recunoască și să evite pericolele.



Indicație

Detalii privind datele de efectuare a pregătirilor, durata și frecvența acestora sunt disponibile la serviciul dvs. local de asistență clienți

Adresa și numărul de telefon sunt disponibile pe internet: <u>http://www.tecan.com/customersupport</u>



# 3. Activități preliminare

## 3.1 Despachetarea și verificarea

## 3.1.1 Verificarea pachetului livrat

Instrumentul livrat include:

- Sursă externă de energie electrică
- Cablu de rețea
- Cablu USB pentru conexiunea la un calculator extern
- Manual de utilizare (IFU), fișiere PDF pe suportul de date
- Stick USB
  - o Software Magellan
  - o Manual de utilizare (IFU), fişiere PDF
  - o Instrumente (de ex. Adobe Reader)





#### PRECAUŢIE

Indicație Pentru a evita pierderea nedorită de date sau atac de viruși/malware, nu înlăturați niciodată protecția la scriere din stick-ul USB.

#### CITITORUL A FOST TESTAT CU CABLUL USB LIVRAT. DACĂ ESTE UTILIZAT UN ALT CABLU USB, NU SE MAI POATE GARANTA FUNCȚIONAREA CORECTĂ A INSTRUMENTULUI.

## 3.1.2 Procedura de despachetare

- 1. Inspectați vizual pachetul dacă prezintă daune, înainte de a-l desface. *Comunicați imediat orice daună constatată.*
- Alegeţi un loc pentru aşezarea instrumentului. Locul trebuie să fie neted, lipsit de vibraţii, ferit de razele soarelui şi fără a conţine praf, solvenţi şi vapori de acizi. Asiguraţi-vă că distanţa dintre instrument şi perete sau orice alt echipament este de cel puţin 5 cm.
- 3. Ridicați instrumentul din ambalajul de carton și așezați-l la locul ales. Aveți grijă, atunci când ridicați instrumentul.
- Inspectaţi vizual instrumentul pentru a observa dacă există piese desprinse, îndoite sau rupte. Raportaţi imediat orice daună.
- Comparaţi numărul de fabricaţie al instrumentului de pe placa de bază a instrumentului cu numărul de fabricaţie de pe avizul de însoţire al mărfii. Raportaţi imediat orice deosebire.
- 6. Verificați accesoriile instrumentului, confruntându-le cu lista pieselor.
- 7. Păstrați ambalajul pentru transportarea ulterioară.

## 3.2 Cerințe de energie

Instrumentul percepe automat tensiunea de alimentare. Prin urmare, nu este necesară modificarea domeniului de tensiune. Verificați specificațiile de tensiune și asigurați-vă că tensiunea de alimentare a instrumentului este corectă conform următoarelor specificații:

#### Tensiune

Instrument de bază cu adaptor AC	100 – 240 V AC, 50/60 Hz
Instrument de bază fără adaptor AC	24 V DC

Dacă tensiunea menționată mai sus nu este disponibilă în țara dumneavoastră, vă rugăm să contactați serviciul local de asistență clienți Tecan.

Conectați instrumentul numai la un sistem de alimentare cu energie electrică, cu legare la pământ de protecție.



#### PRECAUȚIE NU UTILIZAȚI INSTRUMENTUL ÎNTR-UN DOMENIU INCORECT DE TENSIUNE. DACĂ INSTRUMENTUL ESTE PORNIT CU TENSIUNE INCORECTĂ, ACESTA SE VA DETERIORA.



## PRECAUŢIE

NU ÎNLOCUIȚI CABLURILE DE REȚEA DETAȘABILE CU CABLURI CU CARACTERISTICI NECORESPUNZĂTOARE.

## 3.3 Cerințe de mediu înconjurător

Instrumentul trebuie așezat pe o suprafață netedă, plană, lipsită de praf, solvenți și vapori de acizi.

Trebuie evitate vibrațiile și razele directe ale soarelui pentru asigurarea unor rezultate corecte.

Temperatura ambientală:	
Operarea	15 °C - 35 °C (59 °F - 95 °F)
Depozitarea	-30 °C - 60 °C (-30,00 °F - 60,00 °F)
Umiditate relativă:	20 % - 80 % necondensabil la temperatura de funcționare



## 3.4 Cerințe de sistem

	Suportat	Recomandat
PC	PC compatibil Windows cu procesor compatibil Pentium de 1 GHz (Dual Core)	2 GHz (Dual Core)
Sistem de operare	Windows 10 (32 biţi)	
	Windows 10 (64 biţi) Ediţiile: Pro	
Memorie	Windows 10 (32 biţi): 1 GB RAM	2 GB RAM
	Windows 10 (64 biţi): 2 GB RAM	4 GB RAM
Spațiu liber pe hard disk	3 GB	5 GB
Monitor	Super VGA Graphics	
Rezoluție	1024 x 600 și mai mare	1920 x 1080
Adâncime de culoare	256	
Mouse	Mouse Microsoft sau dispozitiv de indicare compatibil	
Comunicație	1 x USB 2.0	2 x USB 2.0 1 x RS232 (serial)
Dispozitive	Windows 10: Dispozitiv grafic DirectX 9 cu WDDM 1.0 sau driver mai recent	
.NET	Microsoft .NET Framework 3.5: În Windows 10 utilizatorului i se va solicita să instaleze versiunea .NET framework necesară (3.5) dacă nu este deja instalată.	
Windows Installer	3.1	
	Dacă această versiune nu este instalată, programul de instalare/actualizare o va instala.	
Microsoft Excel	2007, 2010, 2013, 2016 (32 biţi),	2010 (32 biţi)
	2019 (32 biţi),	2019 (32 biţi)
	Microsoft Excel 365 (32 biţi)	
	Sunt suportate numai edițiile de 32 biți!	
	Edițille Starter NU sunt suportate!	



## 3.5 **PORNIREA** instrumentului

Următoarele proceduri descriu în detaliu etapele necesare, înainte de pornirea instrumentului.



#### PRECAUŢIE

ÎNAINTE DE INSTALAREA ȘI PORNIREA INSTRUMENTULUI, ACESTA TREBUIE LĂSAT CEL PUȚIN TREI ORE ÎNAINTE DE A FI PORNIT, PENTRU A NU EXISTA NICIO POSIBILITATE DE CONDENSARE CARE AR PUTEA CAUZA UN SCURTCIRCUIT.

Atunci când cerințele menționate mai sus au fost respectate, instalarea se execută, utilizând următoarea procedură:

- 1. Conectați instrumentul la calculatorul extern cu cablul USB.
- 2. Asigurați-vă că întrerupătorul general de pe partea stângă a instrumentului este în poziția OFF (OPRIT).
- 3. Introduceți cablul electric în priza de rețea din panoul stâng.
- 4. Porniți instrumentul, utilizând întrerupătorul general de pe panoul stâng.

Instrumentul este pregătit pentru măsurarea microplăcilor după instalarea software-ului.



Indicație Înainte de începerea măsurărilor, asigurați-vă că poziția microplăcii A1 este introdusă corect.

Microplăcile pot fi măsurate numai fără capace.

Închideți capacul suportului plăcii, înainte de a începe o măsurare, pentru a evita influențarea rezultatelor de către lumina ambientală.



Indicație Descărcați întotdeauna microplaca direct de la cititor după finalizarea măsurătorii.



## 3.6 Software

## 3.6.1 Introducere / Domeniu de utilizare

Software-ul de control al instrumentului și de analiză a datelor **Magellan** este furnizat împreună cu instrumentul.

Magellan este **un software universal de control al cititorului şi de analiză a datelor** pentru analizarea datelor generate de testele cu microplăci, folosind dispozitivele de măsură Tecan.

Magellan este disponibil în două versiuni:

- Magellan şi
- Magellan Tracker

Software-ul **Magellan** este destinat analizelor endpoint (punct final), cinetice şi multilabel cu instrumentul INFINITE F50 PLUS în conformitate cu domeniul de utilizare. A se vedea capitolul 2.1 Domeniu de utilizare.

**Magellan Tracker** oferă toată funcționalitatea necesară pentru a se conforma cu Reglementarea FDA 21 CFR partea 11, suplimentar față de funcționalitatea software-ului Magellan.



Indicație

Este important să aveți în vedere că instalarea corespunzătoare a instrumentului și software-ului Magellan nu va asigura conformitatea cu legile și cerințele legale. Principiile corespunzătoare referitoare la procesele și procedurile standard de operare, inclusiv validarea și controlul calității, trebuie de asemenea să fie stabilite.

## 3.6.2 Procedura de instalare

Pentru a instala software-ul, introduceți stick-ul USB în portul USB și procedați astfel:

- Asistentul de instalare a software-ului Magellan trebuie să pornească în mod automat şi să vă ghideze în timpul procesului de instalare. Dacă acest lucru nu are loc, vă rugăm să executaţi fişierul 'E:\Tecan.exe' (E reprezintă litera pentru drive-ul stick-ului USB).
- Selectaţi 'Magellan software' şi în funcţie de versiunea pe care aţi comandato - 'Install Magellan' sau 'Install Magellan Tracker' pentru a începe procedura de instalare şi respectaţi asistentul la instalare.
- 3. Dați clic pe **Instali (Instalare)** pentru a începe procedura de instalare a software-ului.
- 4. Dați clic pe l accept the terms of the license agreement și Next pentru a continua.
- 5. Apare pagina **Customer Information (Informații clienți)**: vă rugăm să introduceți numele utilizatorului și organizația.
- 6. Apare Configuration page (Pagina de configurare): alegeți limba.
- 7. Pagina Use for regulated environments: dați clic pe Next pentru a continua.
- 8. Dați clic pe Install (Instalare) pentru a începe instalarea.
- 9. Dați clic pe **Finish (Terminare)** pentru a încheia instalarea și pentru a alege programul de setare.

#### 3. Activități preliminare



Software-ul poate fi pornit prin meniul Windows **Start**, selectând **Magellan** în grupul de programe **Tecan**.





Indicație Este foarte important ca persoana care instalează software-ul să dețină drepturi de administrator al calculatorului.

Indicație Magellan V7.5 nu poate fi instalat alături de alte versiuni ale softwareului Magellan.

Conform setării iniţiale, toate tipurile de fişiere asociate cu Magellan pentru INFINITE F50 sunt memorate în subdirectorii corespunzători, în următorul director:

- Windows XP:
  - C:\Documents and Settings\All Users\Documents\Tecan\Magellan for F50
- Windows 7, Windows 8, Windows:
   C:\Users\Public\Documents\Tecan\Magellan for F50

## 3.6.3 Calificarea instalării

Verificați instalarea cu succes a software-ului Magellan cu ajutorul programului automat de calificare a instalării:

Start *TecanIQ.exe* din calea inițială de instalare (C:\*Program Files\Tecan\Magellan*) sau din **meniul Windows Start: Start > Programs > Tecan > MagellanIQ.** 

Dați clic pe **Check (Verificare)** pentru a începe calificarea instalării. Toate componentele instalate trebuie să aibă statusul **OK**. Vă rugăm să contactați reprezentantul dumneavoastră local de vânzări Tecan, dacă este raportată orice posibilă problemă.

Pentru a închide programul de calificare a instalării, dați clic pe **Cancel (Anulare)** sau **Exit (leşire)**.



#### Indicatie

Calificarea instalării trebuie repetată ori de câte ori software-ul Magellan este instalat sau actualizat la o versiune mai nouă.



## 3.6.4 Începerea lucrului cu Magellan

Tipul principal al interfeței utilizatorului în software-ul Magellan este asistentul la instalare.

Asistenții Magellan reprezintă module workflow, care sunt ghiduri etapizate pentru executarea unor proceduri complexe.

Ocazional, meniurile sunt disponibile în bara superioară. **Menu (Meniu)** oferă o modalitate convenţională de utilizare a software-ului: item-ul relevant al meniului este selectat din meniurile principale. Toate acţiunile următoare sunt pornite instant sau este afişată o casetă de dialog, unde se pot face selecţii sau intrări suplimentare.

#### Interfaţa utilizatorului - Listă asistenţi

După lansarea software-ului Magellan, apare Wizard List (Listă asistenți) :

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

Fiecare asistent poate fi pornit prin dublu clic sau prin selectarea acestuia și dând

clic pe butonul Next (Continuare)

#### Start Measurement Wizard (Asistent Start măsurare)

Start Measurement wizard (Asistent Start măsurare) include următoarele opțiuni:

- Obtain Raw Data (Obţinerea datelor brute) este utilizată pentru generarea rapidă şi facilă a datelor brute, prin setarea parametrilor necesari de măsură şi prin începerea unei măsurări.
- Use Predefined Method (Utilizare metodă predefinită) este utilizată pentru executarea măsurărilor bazate pe metode definite anterior.
- Start Favorite (Start favorit) este utilizată pentru a selecta una dintre cele mai frecvente metode utilizate din lista de icoane numerotate.

După finalizarea măsurării, se creează un fișier pentru spațiu de lucru.

#### **Evaluate Results Wizard (Asistent Evaluare rezultate)**

**Evaluate Results wizard (Asistent Evaluare rezultate)** este utilizat pentru a vizualiza datele brute și pentru a evalua rezultatele. Parametrii de evaluare pot fi vizualizați și datele pot fi reevaluate.

#### Attach Signature Wizard (Asistent Ataşare semnătură)

Attach Signature wizard (Asistent Ataşare semnătură) este utilizat pentru a semna fişierele de metodă și de spațiu de lucru. Această caracteristică este disponibilă cu Magellan Tracker.

#### Create/Edit a Sample ID List Wizard (Asistentul Creare/Editare listă ID de probe)

Create/Edit a Sample ID list wizard (Asistent Creare/Editare listă ID de probe) este utilizat pentru a crea noi liste ID de probe sau de a edita liste ID de probe, deja existente.

#### Create/Edit a Method Wizard (Asistent Creare/Editare metodă)

Create/Edit a method wizard (Asistentul Creare/Editare metodă) este utilizat pentru a defini sau pentru a edita metode.



Indicație Pentru informații detaliate despre software, vă rugăm să consultați Manualul de utilizare Magellan.

Vă rugăm să aveți în vedere că unele caracteristici descrise în Manualul de utilizare Magellan este posibil să nu fie relevante (disponibile) în Magellan V7.5 în combinație cu INFINITE F50 PLUS. Totuși, toate informațiile necesare sunt descrise în acest Manual de utilizare.



Indicație Vă rugăm să găsiți un exemplu detaliat al unei măsurări ELISA în capitolul 6 Exemplu de aplicație.

Fișierele de metodă Sunrise create cu Magellan V7.x sau mai noi, pot fi deschise cu Magellan V7.5, parametrii de măsurare sunt convertiți în mod automat.

Metodele Sunrise create cu versiuni anterioare trebuie convertite prin folosirea opțiunii 'Convert From' din meniul diverse/fișiere.



## 3.7 Magellan - Editor parametrii de măsură

**Measurement Parameter Editor (Editorul de parametrii de măsură)** este utilizat pentru setare workflow-uri. Fiecare workflow se creează uşor prin selectarea și introducerea etapelor procesuale într-o secvență conform aplicației. Workflow-ul aplicației este apoi vizibil utilizatorului în panoul workflow. Fiecare etapă de proces (element de program) poate fi copiat și lipit (folosind shortcuturile standard Windows Ctrl-C, Ctrl-V sau meniul sensibil la context al mouseului) și mutat în poziția dorită în workflow.

Measurement parameters - infinite F50		
🔶 Lab Ware 🛞	state 1	Selection
Plate     Part of Plate     Measurements	Plate definition: [GRESSR] - Greiner 95 Flat Transporent V Details	Norming selected
Absorbance	PartofPlate A1:H12     2	
Actions (*) Shaking Move Plate	Wavelength         3           Wavelength         405 rm           Reference:         405 rm	
Kinetic <ul></ul>		
Miscellaneous     Image: Comment       User Request     Wait (Timer)       Image: Image: Comment     Image: Comment	Workflow pane	
	(Panou workflow)	Info
(Bară de control)		(Panou info)
Number of plotes: 1 © Help Cancel <<< Back	CHOOSE MEASUREMENT PARAMI	TERS

Measurement Parameter Editor (Editor de parametrii de măsură) constă din următoarele item-uri, care sunt descrise detaliat în capitolele următoare:

- Control Bar (Bară de control)
- Workflow Pane (Panou workflow)
- Info Pane (Panou Info)



## 3.7.1 Control Bar (Bară de control)

**Control bar (Bara de control)** este împărțită în cinci secțiuni. Fiecare secțiune conține elemente de program utilizate pentru a crea un workflow individual.

Puteți crea un workflow, dând dublu clic pe elementul de program selectat sau selectându-l și introducându-l în panoul workflow.

Următoarele elemente de program sunt disponibile, atunci când utilizați INFINITE F50 PLUS:

Lab Ware (Material de laborator)	Plate (Placă) Part of Plate (Parte a plăcii)
Measurements (Măsurări)	Absorbance (Absorbanţă)
Actions (Acţiuni)	Shaking (Agitare) Move Plate (Mutare placă)
Kinetic (Cinetic)	Kinetic Cycle (Ciclu cinetic) Kinetic Condition (Condiție cinetică)
Miscellaneous (Diverse)	Comment (Comentariu) User Request (Solicitare utilizator) Wait (Timer) ((Timp) Aşteptare) Incubation (Incubare)

#### Lab Ware (Material de laborator)

#### Plate (Placă)

Elementul de program **Plate (Placă)** este utilizat pentru a selecta un format de placă din lista drop-down **Plate definition (Definiție placă)**. Dați clic pe **Details (Detalii)**... pentru a vedea informații suplimentare despre placa selectată.

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

#### Part of Plate (Parte a plăcii)

Din setarea iniţială elementul de program **Part of Plate (Parte a plăcii)** este comasat. Atunci când este extins, dând clic, indică o microplacă cu 96 godeuri. Pentru a măsura godeuri individuale, daţi clic pe godeul dorit sau, pentru a măsura un rând de godeuri, selectaţi un cadru în jurul rândului dorit. Daţi clic pe **Details...** (**Detaili)**, pre-vizualizarea plăcii poate fi mărită.

🔹 🔻 Part of Plate	2
1       2       3       4       5       6       7       8       9       10       11       12         A	Details



#### Părți independente ale plăcii

Pot fi selectate părți independente ale plăcii:



Un al doilea rând de godeuri poate fi selectat, apăsând **Control key** de pe tastatură și trasând un cadru peste godeurile, care trebuie selectate.

#### Measurements (Măsurări)

#### Absorbance (Absorbanță)

Elementul de program **Absorbance (Absorbanță)** este utilizat pentru a executa măsurări ale absorbanței. Introduceți sau selectați parametrii respectivi.

Două liste drop-down afişează măsurarea disponibilă și lungimile de undă de referință ale filtrelor, conform filtrelor de absorbanță introduse. Dacă listele dropdown sunt goale, filtrele nu au fost definite.

💧 🔻 Absorb	ance			3
Wavelength-			Label	
Measurement:	492 nm	*	Name: Label1	
Reference:	405 nm	~		

#### Actions (Acţiuni)

#### Shaking (Agitare)

Selectați elementul de program **Shaking (Agitare)** dacă placa trebuie agitată, înainte de măsurare sau între ciclurile cinetice.

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

Introduceți parametrii respectivi:

Duration (Durata)	Introduceți durata procesului de agitare.
Intensity (Intensitate)	Introduceți modul dorit de agitare. Amplitudinea și frecvența sunt afișate, atunci când se alege modul respectiv de agitare.

A se vedea capitolul 4.1.1 Agitare microplacă pentru moduri disponibile de agitare.



#### 3. Activități preliminare

Dând clic pe link-ul <u>Wait a couple of seconds (Aşteptaţi câteva secunde)</u> introduceţi un nou element de program. A se vedea pagina 25 pentru detalii.

#### **Kinetic (Cinetic)**

#### Kinetic Cycle (Ciclu cinetic)

Utilizați elementul de program **Kinetic Cycle (Ciclu cinetic)** pentru a executa mai multe măsurări consecutive, care pot fi executate la anumite intervale.

🎧 🔻 Kinetic Cycle		3
Cycles	C Kinetic Interval	
Number of cycles: 2	Use kinetic interval:	
	⊙ Time: 00:01:00	
0.000	◯ Time: 60000 ≎ ms	
	ns	

Introduceți parametrii respectivi:

Cycles (Cicluri)	Number of cycles (Numărul ciclurilor): Introduceți un număr sau dați clic pe săgețile SUS sau JOS pentru numărul de etape ale măsurării actuale (2 – 1000 cicluri) Duration (Durata): Introduceți durata, format hh:mm:ss.
Kinetic Interval (Interval cinetic)	<b>Use kinetic interval (Utilizați intervalul cinetic)</b> : Introduceți intervalul de timp (hh:mm:ss or ms).

#### Kinetic Condition (Condiție cinetică)

Utilizați elementul de program **Kinetic Condiția cinetică**) pentru a defini ce acțiuni trebuie executate la un anumit ciclu.

🙊 🔻 Kinetic Condition	4
Condition Execute commands at cycle:	

Dacă 2 este introdus pentru **Execute command at cycle (Executare comandă** la ciclul) în cadrul unei măsurări cinetice conţinând, de ex. o etapă de **Shake** (Agitare), agitarea este executată numai la ciclul 2.



Indicație Condițiile cinetice precum Shake (Agitare) trebuie introduse imediat după un element de program Ciclu cinetic, pentru a asigura reproductibilitatea optimă a rezultatului. Utilizatorii sunt sfătuiți să seteze scripturi adecvate înaintea măsurărilor și să folosească același script pentru toate măsurările cinetice similare, pentru a obține rezultate comparabile.

#### **Miscellaneous (Diverse)**

#### **Comment (Comentariu)**

Utilizați elementul de program **Comment (Comentariu)** pentru a introduce o observație sau o afirmație pentru măsurarea actuală în câmpul text.

≢ ▼ Comment	6
Comment:	



#### User Request (Solicitare utilizator)

Elementul de program **User Request (Solicitare utilizator)** informează operatorul instrumentului să execute o acțiune definită în timpul unui workflow, la un anumit moment.

🚯 🔻 User Request	7
Text:	

Dacă de exemplu este utilizat elementul de program **Move Plate (Mutare placă)** pentru a scoate placa în scopul executării unei anumite acțiuni, atunci textul introdus trebuie să informeze operatorul să execute aceste acțiuni. O casetă de dialog indică mesajul și procesul de măsurare se oprește până când se dă clic pe **OK**.

#### Wait (Timer) ((Timer) Aşteptare)

Utilizați elementul de program **Wait (Timer) ((Timer) Aşteptare)** pentru a defini o anumită perioadă de așteptare înainte de următoarea etapă în cadrul executării unui workflow.

Introduceți în câmpul Wait time (Timp de așteptare) timpul necesar.

🮯 🔻 Wait (Timer)		5
Timer       Wait time:     00:01:00       (hh:mm:ss)	Options U Wait for injection I gnore wait at last kinetic cycle	

Introduceți parametrii respectivi:

Timer	Introduceți Timpul de așteptare (hh:mm:ss)
Options (Opţiuni)	Ignorați așteptarea la ultimul ciclu cinetic: Atunci când etapa de programare Wait (Timer) ((Timer) Așteptare) este ultima acțiune în cadrul unei rulări cinetice, timpul de așteptare va fi ignorat în ultimul ciclu.

#### **Incubation (Incubare)**

🮯 ▼ Incubation	4
Timer Incubation time: 00:01:00 😂 (hh:mm:ss)	

Introduceți parametrii adecvați pentru incubare:

Incubation time	Introduceți timpul total (min. 5 s)
(Perioada de	
incubare)	



## 3.7.2 Workflow Pane (Panou workflow)

Workflow pane (Panoul workflow) este fereastra, unde este vizibil scriptul de măsurare și unde sunt definiți și editați parametrii.

Există două modalități de introducere a unui element de program din **Control bar** (Bara de control) în Workflow pane (Panoul workflow):

- Selectaţi un element de program din Control bar (Bara de control); dacă daţi dublu clic pe el, este introdus în Workflow pane (Panoul workflow) imediat după elementul de program anterior.
- Dați clic pe elementul de program din Control bar (Bara de control) și trageți-o în Workflow pane (Panoul workflow) în poziția respectivă.

Elementele de program sunt numerotate conform ordinii lor.

Odată ce un element de program a fost introdus în **Workflow pane (Panoul workflow)**, setările și parametrii pentru acest element pot fi introduși sau editați.

Elementele individuale de program din interiorul **Workflow pane (Panoul workflow)** pot fi comasate pentru a afişa cele mai importante informații sau pot fi extinse pentru a accesa toate funcțiile editabile. Dați clic pe unul dintre triunghiurile de lângă titlul elementului de program, sau , pentru a comuta între două moduri de vizualizare.

Din setarea iniţială, editorul parametrilor de măsură porneşte cu elementul **Plate** (**Placă**), elementul **Part of Plate (Parte a plăcii)** (comasat) și elementul **Absorbance (Absorbanță**) în **Workflow pane (Panoul workflow)**.

Elementele de program selectate actual în cadrul **Workflow pane (Panoul workflow)** sunt afişate cu o linie galbenă la marginea superioară.

Dacă un element de program conține erori sau este invalid în cadrul unui workflow curent, elementul va fi marcat cu un semn de eroare și numărul elementului este evidențiat cu roșu. În **Info pane (Panoul Info)** sunt afișate informații detaliate despre eroare. Dacă workflow-ul conține erori, parametrii de măsură nu pot fi închiși.

#### lerarhia elementelor

lerarhia elementelor în Workflow pane (Panoul workflow) este următoarea:

- 1. Plate (Placă)
- 2. Part of Plate (Range) (Parte a plăcii (rând))

Orice etapă de măsură dorită poate fi introdusă imediat după un element de placă sau rând. Utilizați **Release (Declanşare)** și **Indent (Indentare)** pentru a modifica secvența de executare a componentei individuale de bandă. Selectați un element în **Workflow pane (Panoul workflow)**, dați clic pe butonul drept al mouse-ului și selectați **Release (Declanşare)** sau **Indent (Indentare)**.

Şi alte elemente din **Control bar (Bara de control)** pot fi introduse în ierarhia unui workflow, precum urmează:

Primul element **Range (Rând)** este introdus imediat după elementul **Plate** (**Placă**); apoi pot fi introduse toate elementele următoare **Range (Rând)**.

Etapele cinetice sunt posibile în cadrul unui element **Plate (Placă)** sau **Range** (Rând).

Etapele User Request (Solicitare utilizator), Comment (Comentariu) și Wait (Aşteptare) sunt posibile în cadrul unui element Plate (Placă) sau Range (Rând).



## 3.7.3 Info Pane (Panou Info)

**Info pane (Panoul info)** de pe partea dreaptă a ecranului afişează informații, care sunt relevante pentru elementul de program selectat curent. Sunt afişate orice avertizări și erori.

## 3.8 Magellan - Definirea măsurărilor

Următorul capitol descrie unele exemple pentru a ilustra definirea diferitelor măsurări.

## 3.8.1 Definirea măsurărilor endpoint

Următorul exemplu descrie o Absorbance Endpoint Measurement (Măsurare punct final a absorbanței) în toate godeurile unei microplăci cu 96 de godeuri.

- 1. Selectați o microplacă cu 96 de godeuri din lista drop-down **Plate definițion** (Definiție placă).
- 2. Din setarea iniţială, toate godeurile unei microplăci cu 96 de godeuri sunt alese pentru măsurare.
- 3. Introduceți măsurarea dorită și lungimile de undă de referință.

🔶 🔻 Plate		1
Plate definition:	[COS96R] - Corning 96 Flat Transparent	Details
		Use a part of the plate
🤹 🔻 F	Part of Plate	2
1 2 A C C C C C C C C C C C C C	2 3 4 5 6 7 8 9 10 11 12 Details Details	
4	▼ Absorbance	3
- W	/avelength	
M:	Reference: 620 nm	



## 3.8.2 Definirea măsurărilor multilabel

Următorul exemplu descrie o **Absorbance Multilabel Measurement (Măsurare multilabel a absorbanței)** într-un rând definit al unei microplăci cu 96 godeuri (A1:E7). Trebuie măsurate trei etichete de absorbanță.

- 1. Selectați o microplacă cu 96 de godeuri din lista drop-down **Plate definition** (Definiție placă).
- 2. Din setarea inițială, toate godeurile unei microplăci cu 96 de godeuri sunt alese pentru măsurare.

Dați clic *pentru a extinde elementul Part of Plate (Parte a plăcii)*. Prin urmare, selectați rândul dorit al plăcii (A1:E7).

- 3. Introduceți lungimea de undă dorită a măsurării.
- 4. Introduceți încă 2 elemente **Absorbance (Absorbanță)** și introduceți lungimile de undă ale măsurării.

* Pat of Plate     1     1     2   3   4   5   7   8   9   10 <t< th=""><th></th><th>[COS96it] - Co</th><th>rning 96 Flat Transparent</th><th></th><th></th><th></th><th>Details</th></t<>		[COS96it] - Co	rning 96 Flat Transparent				Details
PatoPlate   1 2 3 4 5 7 8 10 11 12   1 2 3 4 5 6 7 8 10 11 12   1 2 3 4 5 6 7 8 10 11 12   1 2 3 4 5 7 8 10 11 12   1 2 3 4 5 7 8 10 11 12   1 1 2 3 4 5 7 8 10 10   1 1 2 3 4 5 10 10 10 10   1 1 0 0 0 0 0 0 0 0 0 0   1 0 0 0 0 0 0 0 0 0 0 0 0   1 0 0 0 0 0 0 0 0 0 0 0 0 0   1 0 0 0 0 0 0 0 0 0 0 0 0 0 0   1 0 <th></th> <th></th> <th></th> <th></th> <th></th> <th></th> <th> Use a part of the pla</th>							 Use a part of the pla
I 2 3 4 5 6 7 8 9 10 11 12   B I <	🌸 👻	Part of Plate					
Image: Constraint of the state of			7 0 0 40 44 40				
B     F     F     F     F     Kasurement:     Mare:     Label     Wavelength:     Mare:     Label     Wavelength:     Mare:     Label     Wavelength:     Wavelength:	A			Details			
viewelength   Wavelength   Reference:   405 nm     Name:     Labe!     Name:     Labe!     Vavelength   Messurement:   405 nm     Name:   Labe!     Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!   Name:   Labe!     Name:     Labe!     Name:     Labe!     Name:     Labe!     Name:     Labe!     Name:     Labe!     Name:     Labe!     Name:     Labe!     Name:         Name:     Labe!     Name:             Name:	BO		000000				
F   G   H     Vavelength   Name:   Label   Name:   Label   Vavelength   Messurement:   492 nm   Name:   Label   Name:   Label   Vavelength   Name:   Label							
F   G   H     Vavelength   Name:   Label   Name:   Label   Vavelength   Massuement:   405 nm   Vavelength   Name:   Label   Name:   Label   Vavelength   Name:   Label   Name:   Label   Name:   Label   Name:   Label   Name:   Label	E	50000	00000				
S   H     Vavelength   Name:   Label     Name:   Label     Vavelength   Name:   Label   Vavelength   Name:   Label   Vavelength   Name:   Label	F 💽		000000				
Vavelength   Wavelength   Reference:   405 nm   Reference:   405 nm   Name:   Label	G		000000				
Vavelength         Messurement:       405 nm         Reference:       405 nm         Vavelength       Name:         Label       Name:         Vavelength       Name:         Reference:       405 nm         Vavelength       Name:         Reference:       405 nm         Vavelength       Name:         Name:       Label							
Wavelength   Messurement:   405 nm   Reference:   405 nm     Name:   Label							
Wavelength         Reference:         405 nm         Absorbance         Wavelength         Messurement:       492 nm         Reference:       405 nm         Wavelength         Name:       Label         Name:       Label         Wavelength       Name:         Reference:       405 nm         Maxwell       Name:         Label       Name:         Name:       Label		🔻 Absorban	De				
Messurement: 405 nm   Reference: 405 nm     Wavelength   Messurement:   432 nm   Reference:   405 nm     Name:   Label	CV	Wavelength		Label			
Reference:     Wavelength   Messurement:   492 nm   Reference:   405 nm     Name:   Labe!	M	deasurement: 4	05 nm 🔽	Name:	Label1	/	
Wavelength         Messurement:       492 nm         Reference:       405 nm         Wavelength         Varies       Label         Name:       Label		Reference: 4	05 nm 🗸 🗸				
Wavelength       Messurement:       492 nm       Reference:       405 nm       Wavelength       Messurement:       520 nm       Reference:       405 nm							
Wavelength       Label         Measurement:       492 nm         Reference:       405 nm         Wavelength       Label         Measurement:       620 nm         Reference:       405 nm	6	🔻 Absorban	De				
Messurement:     432 nm       Reference:     405 nm       Wavelength       Messurement:     620 nm       Reference:     405 nm	~	Wavelength		Label			
Reference:     405 nm       Wavelength     Label       Reference:     405 nm		vleasurement: 4	92 nm 🗸	Name:	Label2	,	
Wavelength       Reference:       405 nm	M						
Wavelength       Measurement:       E20 nm       Reference:       405 nm	M	Beference: 4	15 pm				
Wavelength Measurement: 620 nm V Reference: 405 nm V		Reference: 4	05 nm 💌				
Measurement. 620 nm V Reference: 405 nm V		Reference: 4	05 nm 🔽				 _
Reference: 405 nm		Reference: 4	05 nm 💌	Label			
		Reference: 4	05 nm 💌	Label	Label3		
		Reference: 4 → Absorban wavelength deasurement: 6 Reference: 4	05 nm 🗸	Label Name:	Label3		
		Reference:     4       ✓     Absorban       Ø     Reference:       4	05 nm 💙 De 20 nm 💌 25 nm 💙	Label-Name:	Label3		

## 3.8.3 Definirea măsurărilor cinetice

Următorul exemplu descrie o măsurare cinetică a unei microplăci cu 96 godeuri.

- 1. Selectați o microplacă cu 96 de godeuri din lista drop-down **Plate definițion** (Definiție placă).
- 2. Introduceți un element de program **Kinetic Cycle (Ciclu cinetic)** între partea plăcii și elementul de absorbanță.
- 3. Cicluri/Numărul ciclurilor: 50
- 4. Interval cinetic (intervalul dintre măsurări): selectați **Use kinetic interval** (**Utilizare interval cinetic**) și introduceți: 2 minute 30 secunde.
- 5. Definiți elementul **Absorbance (Absorbanță)** prin introducerea lungimii de undă dorite a măsurării.

🔷 ▼ Plate	1
Plate definition: [COS96/t] - Corning 96 Flat Transparent	Details Use a part of the plate
🔹 🔻 Part of Plate	2
1       2       3       4       5       6       7       8       9       10       11       12         B <th></th>	
	3
Cycles       So         Number of cycles:       50 \$         Duration       Itime:         0:002:30       (htcmm:ss)         Time:       150000 \$         Time:       150000 \$	
Absorbance	4
Wavelength Measurement: 492 nm V Reference: 405 nm V	



## 3.8.4 Elemente de program Indentare și Declanșare

Decizia de a indenta/declanşa un element de program va modifica workflow-ul unui instrument în timpul măsurărilor.

Acțiunile tuturor elementelor de program cu aceeași indentare sunt executate consecutiv. Singura dependență dintre aceste elemente de program este aceea că următoarea acțiune începe imediat după ce acțiunea anterioară s-a încheiat.

Un element de program care este indentat mai mult decât elementul de program anterior indică dependența dintre cele două elemente de program. Acest lucru înseamnă că parametrii definiți în primul element de program sunt de asemenea activi pentru al doilea element de program (indentat).

Următorul este un exemplu despre definirea unei **Multilabel kinetic (Multilabel cinetice)** cu două **Absorbance labels (Etichete de absorbanță)**. Exemplul indică că două elemente de program **Absorbance (Absorbanță)** depind de elementul de program **Kinetic Cycle (Ciclu cinetic)**, care depinde de elementul de program **Part of Plate (Parte a plăcii)**, care depinde de elementul de program **Plate (Placă)**. Definiți parametrii pentru un exemplu, astfel:

- 1. Plate (Placă): de ex. Greiner 96 plată transparentă
- 2. Ciclu cinetic/Numărul ciclurilor: 5
- 3. Absorbanță/Etichetă lungime de undă 1: 450 nm
- 4. Absorbanță/Etichetă lungime de undă 2: 492 nm

**Workflow pane (Panoul workflow)** apare precum este indicat în screenshot (captură de ecran):

🔷 🔻 Plate		1
Plate definition:	[GRE96it] - Greiner 96 Flat Transparent	Details Use a part of the plate
🄹 🔻	Part of Plate	2
A A B C D C C C C C C C C C C C C C C C C C	2 3 4 5 6 7 8 9 10 11 12 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0	
G	) ▼ Kinetic Cycle	3
	Ouration     Signature         Kinetic Interval         Use kinetic interval	
	🖌 🔻 Absorbance	4
	Wavelength Measurement: 450 nm V Reference: 405 nm V	
	🗼 🔻 Absorbance	5
	Wavelength Measurement: 492 nm M Reference: 405 nm	
		203



Definiția de mai sus rezultă în următorul workflow:

Absorbanța tuturor godeurilor unei microplăci cu 96 godeuri este măsurată prima dată la 450 nm și apoi la 492 nm. Ambele măsurări ale absorbanței sunt executate în 5 cicluri cinetice.

Declanşarea celui de-al doilea element de program Absorbance (Absorbanţă), astfel încât să fie aliniat cu item-ul Kinetic Cycle (Ciclu cinetic), modifică workflow-ul. Selectați al doilea element de program Absorbance (Absorbanţă) și dați clic cu butonul drept al mouse-ului. Selectați Release Strip (Declanşare strip) din meniul sensibil la context. Workflow pane (Panoul workflow) apare precum este indicat în următorul screenshot (captură de ecran):

🔷 ▼ Plate	1
Plate definition: [GRE96it] - Greiner 96 Flat Transparent	Use a part of the plate
🔹 🔻 Part of Plate	2
1       2       3       4       5       6       7       8       9       10       11       12         B       Image: Constraint of the state	
😡 🔻 Kinetic Cycle	3
Cycles     Signature     Kinetic Interval       O Duration     Use kinetic interval	
Absorbance	4
Wavelength Measurement: 450 nm Reference: 405 nm	
🐇 🔻 Absorbance	5
Wavelength Measurement: 492 nm V Reference: 405 nm V	
	203

În acest workflow, o măsurare **cinetică a absorbanţei** cu 5 cicluri este realizată pentru prima absorbanţă la 450 nm; după încheierea acestei bucle, este executată măsurarea **endpoint a absorbanţei** la 492 nm.



## 3.9 Optimizare pentru performanţă maximă

Instrumentul a fost testat în întregime pentru a se asigura că performanţa sa se încadrează în limitele specificate (a se vedea 4.4.2 Specificaţii de măsură pentru detalii).

Instrumentul poate oferi o exactitate maximă, dacă se respectă recomandările indicate mai jos.

### 3.9.1 Amplasarea instrumentului

Instrumentul trebuie amplasat la loc adecvat (a se vedea capitolul 3.3 Cerințe de mediu înconjurător pentru informații detaliate).

### 3.9.2 Procedura de operare

#### Generalități

- Se recomandă respectarea următoarelor proceduri standard de operare pentru analizele utilizate.
- Cea mai bună reproductibilitate se obţine, atunci când lungimea de undă a măsurării corespunde lungimii de undă maxime a absorbanţei a unei soluţii speciale.

Este importantă utilizarea unei lungimi de undă maxime a absorbanţei, în cazul în care curba absorbanţei se află peste o bandă îngustă a lungimii de undă.

Vă rugăm să aveți în vedere că măsurările în pantă a unui vârf al absorbanței va limita acuratețea valorilor OD.

- După ce fiecare microplacă a fost măsurată, vă rugăm să consultați ambalajul setului de testare pentru informații privind procedura de validare.
- Utilizați filtrele recomandate de absorbanță pentru INFINITE F50 PLUS.

#### Microplăci

 Instrumentul poate fi utilizat cu acele tipuri de microplăci care sunt descrise în capitolul 4.4.3 Microplăci. Cele mai bune rezultate sunt obţinute, atunci când sunt utilizate microplăcile cu fund plat. În funcţie de tipul de mircoplacă ce este utilizat, rezultatele de măsură pot diferi.

Aveţi grijă în special atunci când utilizaţi microplăci cu fundul în formă de C, U sau V sau plăci cu stripuri, deoarece este posibil ca rezultatele de măsură să difere faţă de specificaţiile descrise în acest document. Asiguraţi-vă că tipul de microplacă utilizat cu INFINITE F50 PLUS absorbance reader este adecvat pentru aplicaţia respectivă.

- Utilizați numai microplăci perfect curate.
- Nu permiteţi pătrunderea prafului în soluţii sau în microplacă în timpul perioadei de incubare, înainte de măsurare.
   Se recomandă utilizarea unui capac de protecţie la amplasarea unei microplăci în afara instrumentului.
- Inexactitățile în cantitatea de soluție pipetată au un efect semnificativ asupra rezultatelor obținute, atunci când sunt utilizate cantități mici de soluții.
- Forma meniscului soluției poate cauza inexactități în rezultate, în special dacă sunt utilizate cantități mici de soluție.



## 4. Caracteristici instrument

## 4.1 Caracteristici instrument

Următoarele moduri de măsurare a absorbanței sunt disponibile în INFINITE F50 PLUS:

măsurări endpoint, cinetice și multilabel.

## 4.1.1 Agitare microplacă

INFINITE F50 PLUS poate agita microplaca înainte să fie măsurată. Microplaca poate fi agitată între fiecare dintre ciclurile cinetice de măsură.

Utilizați Magellan pentru a seta modul de agitare.



#### PRECAUȚIE ATUNCI CÂND AGITAȚI MICROPLĂCILE, POT APĂREA PIERDERI PRIN SCURGERE, DACĂ GODEURILE SUNT UMPLUTE PREA MULT.

Moduri de agitare pentru INFINITE F50 PLUS:

Mod agitare	Amplitudine agitare	Frecvență agitare
HIGH (ÎNALT)	2,8 mm	12,3 Hz
NORMAL (NORMAL)	4,4 mm	9,2 Hz
LOW (SCĂZUT)	4,4 mm	7,8 Hz
WIDE (AMPLU)	14,1 mm	2,0 Hz

## 4.2 Descriere instrument

Figura de mai jos indică componentele instrumentului.





#### 4. Caracteristici instrument

LED-ul de stare oferă informații despre starea instrumentului:

- Luminare intermitentă pe fond verde: instrumentul nu este conectat la Magellan
- Verde: instrumentul este conectat și pregătit pentru măsurare
- Roşu: măsurarea este în curs

Pe panoul stâng al instrumentului, se pot găsi portul USB, întrerupătorul principal și priza de rețea.



Întrerupător principal

Placa cu nume este ataşată pe partea inferioară a instrumentului.

Exemplu de plăcuță de identificare



Conținutul plăcuței de identificare (de exemplu, numele modelului și numărul articolului) poate varia în funcție de modelul specific.

Pentru o prezentare generală a diferitelor instrumente pentru care este valabil prezentul manual de utilizare consultați Declarația de conformitate ("Declaration of Conformity") de pe ultima pagină a acestui document.



## 4.3 Descriere centrifugă cu filtre

Centrifuga standard cu filtre INFINITE F50 PLUS este furnizată cu patru filtre de interferență cu benzi înguste care au o lungime fixă de undă (405, 450, 620 şi 492 nm). Este posibilă echiparea centrifugei cu până la 8 filtre. Pentru filtrele disponibile ca accesorii, vă rugăm să contactați reprezentantul dumneavoastră local de vânzări Tecan.

Filtrele centrifugei standard cu filtre sunt montate precum este indicat mai jos:

Poziție filtru	Lungime de undă a filtrului
1	405 nm
2	450 nm
3	620 nm
4	492 nm
5 - 8	poziții goale de filtru



Atunci când o lungime de undă este selectată pentru măsurare, filtrul specific este adus în dreptul fasciculului de lumină, mutând centrifuga cu filtre în poziția corespunzătoare.



Indicație Pentru informații suplimentare despre definiția unui nou filtru, a se vedea 7.5.2 Definirea filtrelor.

## 4.4 Specificații instrument

Tabelele de mai jos listează specificațiile pentru INFINITE F50 PLUS absorbance reader.

## 4.4.1 Specificații generale

PARAMETRII	CARACTERISTICI		
Consum de putere Alimentare externă de la reţea	Alimentare de la reţea: Instrument de bază cu adaptor AC: 100-240 V AC, 50/60 Hz, max. 1.2 A (auto sensing, supratensiune categoria II) Instrument de bază fără adaptor AC: 24 V DC (supratensiune categoria I)		
Consum INFINITE F50 PLUS	Mod de aşteptare: aprox. 12 VA Mod operațional: max. 30 VA		
Dimensiuni exterioare	Lăţime 34,7 cm(13,66 inch) Adâncime: 18,9 cm( 7,44 inch) Înălţime: 13,4 cm( 5,28 inch)		
Greutate	2.6 kg (inclusiv alimentarea de la reţea)		
Temperatura ambientală:			
Operarea	15 °C - 35 °C (59 °F - 95 °F)		
Depozitarea	-30 °C - 60 °C (-30,00 °F - 60,00 °F)		
Umiditate relativă	20 % - 80 %		
Grad de poluare	2		
Metodă de eliminare	Deşeuri contaminate		
Mediu înconjurător	A se vedea capitolul 3.3 Cerințe de mediu înconjurător pentru informații suplimentare.		


# 4.4.2 Specificații de măsură

PARAMETRII	CARACTERISTICI
Timp de măsură: o singură lungime de undă lungime de undă dublă	< 15 secunde < 20 secunde
Interval lungime de undă: Standard	400 - 750 nm
Interval lungime de undă: 400 - 750 nm	0 - 4.000 OD
Rezoluție:	0.0001 OD
Exactitate: 450, 492 nm 0.000 - 2.000 OD 2.000 - 3.000 OD	≤ (1.0 % + 0.010 OD)* ≤ (1.5 % + 0.010 OD)*
Precizie: 450, 492 nm 0.000 - 2.000 OD 2.000 - 3.000 OD	≤ (0.5 % + 0.005 OD)* ≤ (1.0 % + 0.005 OD)*
Linearitate: 450, 492 nm 0.000 - 2.000 OD 2.000 - 3.000 OD	≤ 1 % ≤ 1.5 %
Selectare lungime de undă: Filtru standard	Filtre de interferență cu bandă îngustă. Într-o centrifugă cu filtre pot fi montate până la 8 filtre.
Exactitatea lungimii de undă a filtrului:	Lungime de undă centrală ± 2 nm
Lăţime bandă de filtru: La 50 % transmisie	10 ± 2 nm
Sursă de lumină:	LED
Interfață calculator:	USB

Toate dispozitivele conectate trebuie aprobate și listate ca pentru standardele de siguranță și locale echivalente ale Echipamentelor de Tehnologie a Informațiilor IEC 60950-1.

 $^{\ast}$  mai bună decât sau egală cu x % din valoare de măsurare plus valoarea OD corespunzătoare

# 4.4.3 Microplăci

Toate microplăcile cu 96 godeuri cu fundul transparent (plat, sub formă de C, U şi V; inclusiv microplăci cu stripuri) care sunt în conformitate cu următoarele standarde, pot fi utilizate cu INFINITE F50 PLUS absorbance reader:

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



#### PRECAUȚIE UTILIZAȚI NUMAI MICROPLĂCI FĂRĂ CAPACE ȘI NU UTILIZAȚI MICROPLĂCI MAI MARI DE 15.2 MM.



#### Manevrarea microplăcii

Introduceți sau scoateți microplaca numai atunci când suportul plăcii este complet scos (vezi figura de mai jos) și motorul suportului plăcii nu este activ. Nu desfaceți capacul carcasei atunci când LED-ul de stare este roșu.



## AVERTIZARE PURTAŢI MEREU MĂNUŞI RECICLABILE ŞI ÎMBRĂCĂMINTE DE PROTECŢIE ATUNCI CÂND MANEVRAŢI MICROPLACA.



Poziția godeului A1

Suport al plăcii – complet scos

# 4.5 Accesorii instrument

Lista de mai jos conţine accesoriile opţionale disponibile pentru INFINITE F50 PLUS, care pot fi comandate suplimentar:

- Filtre suplimentare
- Unealtă de asamblare a filtrelor
- Placă MultiCheck<sup>™</sup> pentru familia INFINITE F50

Pentru informații suplimentare și disponibilitatea în țara dumneavoastră, vă rugăm să contactați reprezentantul dumneavoastră local de vânzări Tecan.



# 5. Control al calității

# 5.1 Introducere



#### PRECAUȚIE DACĂ PERFORMANȚA ANALITICĂ A DISPOZITIVULUI INFINITE F50 PLUS ESTE PUSĂ ORICÂND LA ÎNDOIALĂ, RESPECTAȚI INSTRUCȚIUNILE DATE PENTRU CONTROLUL CALITĂȚII SAU CONTACTAȚI SERVICIUL LOCAL DE ASISTENȚĂ CLIENȚI TECAN.

Acest capitol oferă informații despre procedura de auto-verificare pentru instrument și instrucțiuni despre cât de facil este să verifici calitatea operațională.

# 5.2 Procedura de auto-verificare

În timpul conectării a INFINITE F50 PLUS la software-ul de control al cititorului Magellan, se verifică motoarele și senzorii, și se inițializează suportul plăcilor și centrifuga cu filtre.

Înainte de fiecare măsurare, se execută o procedură de calibrare de autoverificare, pentru a se asigura că instrumentul funcționează corect și pentru a calibra sistemul optic.

# 5.3 Calificare operațională (OQ)

Următoarele teste pot fi executate pentru a se asigura că instrumentul funcționează în mod corect și că se obțin rezultate exacte.

Reproductibilitatea și exactitatea instrumentului pot varia în funcție de tipul de soluție și de mircoplaca utilizată.

Pentru a elimina acest efect, instrumentele sunt testate în fabrică cu o placă de calibrare, care îndepărtează influența soluției și orice variație cauzată de așezarea microplăcii, atunci când este măsurată.

# 5.3.1 Test MultiCheck

Testul Multicheck asigură o verificare automată a performanţei cititorului inclusiv a exactităţii, linearităţii, preciziei şi trasabilităţii conform standardelor furnizate de NIST.

# 5.3.2 Testarea microplăcii

Dacă densitățile optice ale godeurilor din microplacă nu sunt consistente, rezultatele obținute cu acest tip de mircoplacă va fi influențat.

Inconsistența poate fi verificată citind o microplacă goală.

Valorile OD obținute din măsurarea unei microplăci goale ar trebui să se găsească într-un domeniu îngust. De exemplu: ± 0.010 OD.

Dacă valorile OD nu se încadrează în acest domeniu, acest tip de microplacă nu trebuie utilizată.

Prin folosirea măsurărilor duale ale lungimii de undă, influența diferenței în valorile OD ale microplăcii este îndepărtată sau redusă la un nivel, care se încadrează în limitele acceptabile.



# 5.3.3 Precizia instrumentului cu probele de lichid

Această procedură poate fi utilizată pentru a verifica reproductibilitatea măsurărilor. Se recomandă utilizarea unei microplăci cu fund plat.

Umpleți o nouă microplacă cu o soluție Orange G proaspăt preparată; utilizați diferite diluări ale soluției în fiecare godeu, astfel încât să se obțină o gamă de densități optice. Asigurați-vă că fiecare godeu conține cel puțin 200 µl. Seriile de diluare trebuie să se încadreze în domeniul 0.1 - 3.0 OD. Pentru a obține aproximativ 3 OD, se recomandă utilizare a 125 mg.l<sup>-1</sup> de soluție Orange G (Sigma, Nr. cat. O7252).

Definiți o funcționare de probă, utilizând un filtru de 492 nm și apoi măsurați microplaca de cel puțin trei ori.

Calculați pentru fiecare godeu:

- media valorii OD
- deviația standard

## Exemplu

#### Citiri între 0.000 - 2.000 OD

Deviația standard a fiecărui godeu trebuie să se încadreze (0.5 % + 0.005 OD).

Calcularea deviației maxim admise, utilizând 1.000 OD, ca valoare OD medie:

1.000 \* 0.5 % + 0.005 = 0.010 OD

#### Citiri între 2.001 - 3.000 OD

Deviația standard a fiecărui godeu trebuie să se încadreze (1,0 % + 0.005 OD).

Calcularea deviației maxim admise, utilizând 2,400 OD, ca valoare OD medie:

2,400 \* 1,0 % + 0.005 = 0,029 OD

#### Citiri peste 3.000 OD

Citirile care depăşesc 3.000 OD sunt utilizate numai ca o indicație și precizia nu poate fi garantată.

# 5.3.4 Linearitatea instrumentului cu probe de lichid

Linearitatea instrumentului și aplicația la lungimea de undă utilizată pot fi verificate utilizând o serie de diluări ale unei soluții.

Rezultatul depinde de puritatea colorantului utilizat și de meniscul lichidului din godeuri.

Ca referință, poate fi utilizată o serie de diluări ale soluției Orange G pentru măsurări la 492 nm.

Seriile de diluare trebuie să se încadreze în domeniul 0.1 - 3.0 OD. Pentru a obține aproximativ 3 OD, se recomandă utilizare a 125 mg.l<sup>-1</sup> de soluție Orange G (Sigma, Nr. cat. O7252).

Pentru alte lungimi de undă, trebuie utilizate diferite soluții.

200 µl din fiecare diluare sunt pipetați în microplacă, trebuie utilizat un minim de cel puțin două probe pentru fiecare diluare pentru a reduce erorile cauzate prin pipetare.

Microplaca este apoi măsurată și se trasează o regresie liniară de OD contra concentrației din media valorilor OD măsurate.

Determinați valoarea reziduală neevaluată la pătrat R<sup>2</sup> a liniei de regresie.



Valorile reziduale tipice la pătrat pentru o aplicație standard sunt egale sau mai mari decât  $R^2 = 0.998$ .



Indicație Datele pot varia din cauza inexactității de pipetare.



# 6. Exemplu de aplicație

# 6.1 Introducere

**Fişierele exemplu** Magellan oferă metode Magellan și spații de lucru pentru prezentarea software-ului și pentru a-i facilita utilizatorului munca cu acesta. Fișierele exemplu pentru o analiză cantitativă și calitativă ELISA sunt instalate în mod automat după instalarea software-ului Magellan.

# 6.2 Exemplu detaliat: Analiză cantitativă ELISA

Un exemplu detaliat (test cantitativ) despre modul de crearea a unei metode în Magellan este prezentat în acest capitol. Urmând instrucțiunile, veți învăţa cum să definiți evaluările dintr-o descriere a setului de testare în Magellan.



Indicație Fișierele exemplu apar în mod automat în Lista metodelor din Magellan. Pentru Magellan Tracker, aceste fișiere sunt disponibile din setarea inițială și trebuie convertite.

# 6.2.1 Descrierea setului de testare

În descrierea setului de testare al producătorului a unei ELISA cantitative de detectare a anticorpilor IgM, se găsesc următoarele instrucțiuni: Plate Layout (Dispunere placă)

	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										

BLK = Blank, NC = Negative control, C1 - C6 = Aparate de etalonare (Standarde), S1 - S... = Probe



#### 6. Exemplu de aplicație

#### Măsurare și evaluare

Citiți placa la o lungime de undă de 492 nm, referință la 620 nm.

Cititor valoare blank/placă la godeul A1.

Concentrațiile aparatelor de calibrare (standarde):

Aparat de calibrare 1	5 UA/ml
Aparat de calibrare 2	10 UA/ml
Aparat de calibrare 3	20 UA/ml
Aparat de calibrare 4	40 UA/ml
Aparat de calibrare 5	80 UA/ml
Aparat de calibrare 6	160 UA/ml

După corecția valorii blank, densitățile optice (OD 492 - OD 620) sunt măsurate în comparație cu concentrația. Linia de regresie, care trece prin aceste puncte este curba standard.

Interpretarea rezultatelor testului:

IgM < 18 UA/ml	Negative (Negativ)
18 UA/mI ≤ IgM < 22 UA/mI	Intermediate (Intermediar)
IgM ≥ 22 UA/mI	Positive (Pozitiv)

Concentrația IgM calculată a ambelor controale negative trebuie să se afle sub 8 UA/ml.

#### **Prelucrarea datelor**

În urmă măsurării, fişierul de date (spaţiu de lucru) este stocat în mod automat şi este creat un raport care conţine parametrii de măsură, amplasarea plăcuţei, valorile blanked, curbă standard, concentraţii IgM, definiţia valorii critice, rezultatele calitative ale probelor şi validărilor.

În plus, amplasarea și rezultatele calitative sunt stocate ca fișier ASCII.



# 6.2.2 Crearea unei metode

În caseta de dialog Wizard List (Listă asistenți), selectați Create/edit a method (Creare/Editare metodă) și dați clic pe OK. Dați clic pe Continue (continuare) de pe pagina Welcome (de întâmpinare) a Create/edit a method wizard (asistentului Creare/Editare metodă) și apare caseta de dialog Select a file (selectați un fișier). Selectați New (Nou).

Create/Edit a Method			
New ⊙ Open Control internot	Show:	Files from this instrument  Remarks	Print Preview Status
Help Cancel <<< Back		MAKE YOUR SELE	CTION
		© 2009	Tecan

## Parametrii de măsură

Dați clic pe Make Your Selection (Selectați) și apare pagina Measurement parameters (Parametrii de măsură).

Measurement parameters - infinite F50			
🔷 Lab Ware 🔹	🔷 ▼ Plate	1 Selection	n 🗠
+ Plate	Plate definition: [GRE36it] - Greiner 96 Flat Transparent	V Details	elected
Part or Hate		Use a part of the plate	
Measurements 🔹	🌸 🕨 Part of Plate A1:H12	2	
Absorbance	👃 🔻 Absorbance	3	
Actions 🛞	Wavelength		
🔀 Shaking	Measurement: 405 nm Vame: Label1 V		
Move Plate	Reference: 405 nm		
😱 Kinetic 🏾 🛞			
Kinetic Cycle     Kinetic Cycle			
> Intele contailon			
Miscellaneous 🛞			
Comment User Request			
🍪 Wait (Timer)			
Incubation			
Number of plates: 1			
Help	ſ	HOOSE MEASUDEMENT DADAMETEDS	
Cancel <<< Back		HOOSE MEASUREMENT PARAMETERS	
© 2009 Teran			



# 6. Exemplu de aplicație

Pe stripul **Wavelength (Lungime de undă)** selectați 492 nm ca Measurement wavelength (Lungime de undă a măsurării) și 620 nm ca Reference wavelength (Lungime de undă de referință).

🖕 🔻 Absorb	ance		3
-Wavelength-		Label	
Measurement	492 nm 💌	Name: Label1	
Reference:	620 nm 👻		

Continuați Wizard, dând clic pe **Choose measurement parameters (Alege parametrii de măsură)** și se afișează fereastra **Plate layout (Dispunere plăcuță)**.

Create/Edit a Method														×
Undo Redo Select all unused	10% 1	00% Zoo	om											
Method layout S Plate layout														
cdit method		1	2	3	4	5	6	7	8	9	10	11	12	
	A											Identifiers	Exp. group:	
	В											ST PC	ID-Num 4 1	
	C											LPC HPC BF	Replicates H ③ Fix numb	er.
	D											BF Def. identi		
	E											selection	selection	
	F													
	G													CAN Austri
	H													TEC
Help Cancel <<< Back												© 2009	NEXT	



# **Design Layout**

Definiți dispunerea plăcii, utilizând căsuța de dialog **Well Assignment (Alocare godeu)** de pe partea dreaptă a ecranului.

We	ll Assig	nm.	
r I	dentifier	s	Exp. group:
	SM	^	1
	BL		-ID-Num.
	ST		≠ 4
	PC		
	NC		
	LPC		Replicates
	HPC		≠ 4
	BF		• Fix number:
	RF	~	1
	Def. ide	ntif.	O All
	Fill selecti	on	Delete selection

În caseta de grup Identifiers (Identificatori), selectați BL (Blank).

În caseta de grup Experimental rămâne numărul 1.

În caseta de grup **Replicates (Reproduceri)**, **All (Toate)** este selectat în mod automat.

Dați clic pe godeul A1, care este marcat apoi cu o margine roșie.

Dați clic pe **Fill selection (Selectare umplere)** și godeul este etichetat cu tipul selectat de identificator.



Indicație Un singur godeu poate fi de asemenea umplut, dând dublu clic pe el.

Acum alegeți următoarele setări în caseta de dialog Well Assignment (Alocare godeu):

În caseta de grup **Identifiers (Identificatori)**, selectați **NC** (Negative Control/Control negativ).

În caseta de grup Experimental rămâne numărul 1.

All (Toate) din caseta de grup Replicates (Reproduceri) este selectat în mod automat.

Începând la godeul **B1** dați clic și trageți cu mouse-ul la **C1**. Godeurile **B1** până la **C1** sunt marcate apoi cu o margine roșie.

Dați clic pe **Fill selection (Selectare umplere)** și godeurile sunt etichetate cu tipul selectat de identificator.

Aparatele de calibrare (standarde) trebuie alocate godeurilor **D1** până la **G2**. Alegeți următoarele setări în caseta de dialog **Well Assignment (Alocare godeu)**:

În caseta de grup Identifiers (Identificatori), selectați ST (Standard).





În caseta de grup Experimental rămâne numărul 1.

În caseta de grup **Replicates (Reproduceri)**, alegeți între **Fix number (Stabilire număr)** și **All (Toate)**.

Fix number (Stabilire număr):

Activat numai pentru standarde și probe, unde pot fi utilizate ID-uri. Dacă acest buton **Fix number (Stabilire număr)** este un număr activ, poate fi introdus în câmpul text corespunzător. Acest număr definește cât de multe reproduceri sunt avute în vedere pentru această metodă. În godeurile selectate, este creat numărul introdus de reproduceri pentru fiecare ID. Prin urmare, numărul godeurilor selectate trebuie să fie un multiplu al numărului de reproduceri introdus.

#### All (Toate):

Toate godeurile selectate sunt definite ca reproduceri. Dacă este ales un număr existent de ID pentru probe și standarde, godeurile selectate sunt apoi adăugate ca reproduceri la reproducerile existente. Cu toate celelalte tipuri de identificatori, godeurile selectate sunt adăugate ca reproduceri la reproducerile existente.

Două butoane săgeți definesc direcția reproducerii și secvența numărului ID (orizontal sau vertical).

În acest exemplu, selectați Fix number (Stabilire număr) și 2.

În caseta **ID-Number (Număr ID)** și în caseta de grup **Replicates (Reproduceri)** selectați **vertical arrows (săgeți verticale)**.

Selectați apoi godeurile D1 - G2 și dați clic pe **Fill selection (Selectare umplere)**.



Indicație

Selectați godeurile astfel: Începând la godeul D1, dați clic și selectați cu mouse-ul godeurile necesare până la H1. Apoi țineți apăsată tasta control key (Ctrl) și selectați cu mouse-ul godeurile necesare începând de la A2 până la G2.



Create/Edit a Method															×
Undo Redo Select all unus	sed (	10%	100% Zo	om											
Method layout O	Create/														
Conc, Dil, Refvalues	edit method		1	2	3	4	5	6	7	8	9	10	11 Well Assign	12	
Concentrations  Standard curve		A	BL1 1/1	ST1_3 2/2									Identifiers SM	Exp. grou	
Cutoff definition QC Validation		B	NC1 1/2	ST1_4 1/2									ST PC	-ID-Num 	<b>1</b>
Data handling (©) Data export Printed report		C	NC1 2/2	ST1_4 2/2									LPC	Replicate	
Automated data handing     Miscellaneous     User prompts		D	ST1_1 1/2	ST1_5 1/2									RF Def. ident	it. All	nber:
Number format Method notes		E	ST1_1 2/2	ST1_5 2/2									Fill	Delete	
		F	ST1_2 1/2	ST1_6 1/2											
		G	ST1_2 2/2	ST1_6 2/2											N Austria
		H	ST1_3 1/2												TEC/
Help Cancel <<< Back														NEXT	

Plate Layout (Dispunere placă) apare astfel:

Dați clic pe **Select all unused (Selectare toate neutilizate)** din bara de instrumente pentru a selecta toate godeurile goale de pe placă. Apoi, mențineți apăsată tasta control key (Ctrl) și dați clic pe godeul **H12**, astfel încât să rămână blank și nemarcat.

În caseta de dialog **Well Assignment (Alocare godeu)** selectați **SM (Probă)** de la *Identifiers (Identificatori)*.

În caseta de grup Experimental rămâne numărul 1.

În caseta de grup **Replicates (Reproduceri)**, alegeți **Fix number (Stabilire număr)** și **2**.

În caseta ID-Number (Număr ID) lăsați 1 și în caseta de grup Replicates (Reproduceri) selectați vertical arrows (săgeți verticale). Dați clic apoi pe Fill selection (Selectare umplere). Procedura layout definition (definire dispunere) este completă.

### Transformări

În bara de control de pe partea stângă a ferestrei, selectați următoarea opțiune, Add new transformation... (Adăugare transformare nouă...) din item-ul Transformed data (Date transformate) pentru a defini reducerea valorii blank.

Apare o casetă de dialog, care vă întreabă dacă doriți să definiți reducerea valorii blank. Dați clic pe **Yes (Da)**. Apare următoarea fereastră:



## 6. Exemplu de aplicație

Create/Edit a Method															
Undo Redo	_	10%		om											
<u>ې</u>	)	Input data	Difference data				~							Constants	Options
Method layout 🛞	Creat	f	x <mark>xBL1</mark>							~	Available dat	a (multiple data	sets) 🔻	Functions&Co	nstants 🗨
Conc.+, Dil.+, Refvalues	e/edi														
http://www.transformed.data 🔗	tmet														
Blank reduction Add new transformation	pod	0	) 1	2	3	4	5	6	7	8	9	10	11	12	
Soncentrations			BL1	ST1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37	
Standard curve			1/1	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Evaluate data 🛞			NC1	ST1 4	SM1 2	SM1 6	SM1 10	SM1_1/	SM1 18	SM1 22	SM1 26	SM1 30	SM1 34	SM1 38	
QC Validation		E	3 1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
🏄 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	
Data export		C	NC1 2/2	ST1_4 2/2	SM1_2 2/2	SM1_6 2/2	SM1_10 2/2	SM1_14 2/2	SM1_18 2/2	SM1_22 2/2	SM1_26 2/2	SM1_30 2/2	SM1_34 2/2	SM1_38 2/2	
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	
🛞 Miscellaneous 🛞			ST1_1	ST1_5	SM1_3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
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	
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	
			2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
			ST1 2	ST1 6	SM1 4	SM1 8	SM1 12	SM1 16	SM1 20	SM1 24	SM1_28	SM1 32	SM1 36	SM1 40	
		6	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	m.
		G	2/2	2/2	5M1_4 2/2	SM1_8 2/2	5M1_12 2/2	SM1_16 2/2	SM1_20 2/2	SM1_24 2/2	SM1_28 2/2	SM1_32 2/2	SM1_36 2/2	2/2	Austr
			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
		l n	ST1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37		Ĕ
			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1		
Help														NEYT	
Cancel (<<< Back															
													(D) 200	19 Tecan	

Difference Data (Date diferențiale) este selectat în mod automat în caseta Input data (Date de intrare). Dacă ați confirmat anterior definiția reducerii valorii blank, software-ul îl numește în mod automat Blank reduction (Reducere valoare blank) (vezi datele transformate din bara de control).

În caseta **Formula (Formulă)** apare în mod automat **x-BL1** pentru această reducere a valorii blank, unde x se referă la valoarea curentă a datelor de intrare dintr-un godeu și BL1 este valoarea medie a godeurilor blank ale grupului experimental 1.

Pentru detalii și explicații suplimentare referitor la definirea și alocarea transformărilor, consultați Manualul de utilizare Magellan.

În fiecare godeu, apar următoarele informații (exemplu godeu A5):

SM1_9	Probă, număr grup experimental 1, număr ID probă 9.
2/2	Numărul reproducerii este 2, numărul total al reproducerilor este 2.
x-BL1 sau 1	Transformare alocată x-BL1 (atunci când este selectată Transformation/Transformare) sau Valoare a Factorului de Diluție 1 (atunci când este selectat valori Conc., Dil,. Ref.).



# Definirea valorilor de Concentrație / Diluție / Referință

În bara de control, selectați **Conc., Dil., Ref.-value**s (Valori Conc., Dil., Ref.) din item-ul **Method layout (Configurare metodă)** pentru a defini valorile respective așa cum sunt descrise în setul de testare.

Aparat de calibrare 1	5 UA/ml
Aparat de calibrare 2	10 UA/ml
Aparat de calibrare 3	20 UA/ml
Aparat de calibrare 4	40 UA/ml
Aparat de calibrare 5	80 UA/ml
Aparat de calibrare 6	160 UA/ml

Asigurați-vă că ST este selectat în lista Select Identifier (Selectare identificator).

În lista **Identifier (Identificator)** apare lista standardelor din Grupa Exp. 1. În caseta corespunzătoare **Concentration (Concentrație)** a tipului **ST1\_1**, numărul **5** și în caseta **Unit (Unitate)**, tipul UA/ml. În caseta corespunzătoare **Concentration (Concentrație)** a tipului **ST1\_2** numărul **10**. Unitate trebuie să fie definită numai o singură dată și este valabilă pentru toate standardele. Introduceți valorile pentru ST1\_3 - ST1\_6 în același mod.

Este afișat ecranul care indică dispunerea plăcii și concentrația:





# Standard Curve (Curbă standard)

În bara de control, dați clic pe **Standard curve (Curbă standard)** din item-ul **Concentrations (Concentrații)** pentru a defini curba standard corespunzătoare.

Următoarea se află în descrierea setului de testare:

După corecția valorii blank, densitățile optice (OD 492 - OD 620) sunt măsurate în comparație cu concentrația. Linia de regresie, care trece prin aceste puncte este curba standard.

Pe bara **Data (Date)**, selectați **Blank reduction (Reducere valoare blank)** ca date de intrare.

Data	Analysis type	Intercepts Axis Graph
	1	
	input data:	Blank reduction
	⊙ Standards fro	/m Layout
	<ul> <li>Standards from</li> </ul>	m ext. file: Select
	⊖ Standards fro	m experimental group:
	🔿 No Standard	curve graph
	Additional c	oncentrations

Pe bara **Analysis type (Tip analiză)**, selectați Linear regression (Regresie lineară).

Point to point     Inear regression     Non-linear regression     Lubic spline     Akima	Data scaling: Lin(x)Lin(y)
<ul> <li>Polygomial</li> <li>Eour parameters</li> <li>Eour parameters Marquardt</li> <li>Five parameters</li> <li>LogitLog</li> </ul>	More
Include (0.0)	



Pe bara **Axis (Ax),** definiți etichetarea și scalarea axului, precum este indicat mai jos:

l abel:	Concentration [LIA/m]	
0.1		
Lolor:		Log. Scaling
💿 Auto	select range	
🔘 Rang	e	Min.: Max.:
🔽 Grid		Color: Style:
r'-axis		
Label:	Blank reduction	
Color:	<b>•</b>	Log. Scaling
💿 Auto	select range	
🔘 Rang	e	Min.: Max.:

Pe bara Graph (Grafic), definiți titlul, curbele, fontul grafic și afișajul grafic.

Data	Analysis type Intercepts Axis Graph							
	Tile							
	Label: IgM-ELISA							
	Curves							
	Labet Gro 1							
	Color							
	Symbol							
	Size: Medium 🕑							
	Font Display							
	⊙ Small ✓ Legend ✓ Intercepts							
	○ Medium     Image       ○ Large     Image							



#### 6. Exemplu de aplicație

#### **Definire valori critice**

În bara de control selectați **Cutoff definition (Definire valoare critică)** din itemul **Evaluate data (Evaluare date)** pentru a defini limitele evaluării calitative.

Descrierea setului de testare conține următoarele instrucțiuni:

Interpretarea rezultatelor testului:

IgM < 18 UA/mI	Negative (Negativ)
18 UA/ml ≤ IgM < 22 UA/ml	Intermediate (Intermediar)
IgM ≥ 22 UA/mI	Positive (Pozitiv)

Utilizați următoarea procedură pentru a defini valorile critice adecvate:

În caseta Date de intrare, selectați Mean conc. (UA/ml) (Conc. medie).

Tabelul **Cutoffs (Valori critice)** reprezintă o scală care indică capătul superior și inferior pentru **Limits (Limite)** și **Labels (Etichete)**. În **Limits (Limite)**, introduceți 22 ca prima limită (superioară) și 18 ca a doua limită (inferioară).

În Labels (Etichete), introduceți interpretarea testului (Positive (Pozitiv), Intermediate (Intermediar) și Negative (Negativ)) în casete individuale. Utilizați paleta de culori drop-down pentru alocarea unei culori:

- Positive (Pozitiv) Roşu
- Intermediate (Intermediar) Albastru
- Negative (Negativ) Verde

Ecranul conține următoarele:

▼ positive		
 <ul> <li>intermedial</li> </ul>	10	 
<ul> <li>negative</li> </ul>		
<u> </u>		
<u> </u>		
	Low	
	ļĂ	

Dați clic pe **Cutoff results selection (Selectare rezultate valori critice)** pentru a selecta tipurile de identificatori pentru care trebuie indicate rezultatele valorilor critice.



# Definire QC Validations (Validări QC)

În bara de control, dați clic pe **QC Validations (Validări QC)** din item-ul **Evaluate data (Date evaluare)**. Trebuie definite criteriile de validare pentru test, astfel încât să fie garantată valabilitatea rezultatelor testului.

În acest exemplu, trebuie îndeplinită următoarea cerință:

Concentrația IgM calculată a ambelor controale negative trebuie să se afle sub 8 UA/ml.

În caseta Input, selectați Single conc. (UA/ml) (Conc. unică).

Scrieți în primul rând NC1\_1<8



Indicaţie NC1\_1 înseamnă Control negativ al grupului experimental 1, reproducere 1.

Scrieți în al doilea rând NC1\_2<8

Caseta de dialog QC Validations (Validări QC) este afișată acum astfel:

the second second second second	0.1								
input uata.	Single cor	nc. (UA/I	mij			~		Validation group:	
Validation C	onditions	:							<u>^</u>
NC1_1<8									
NC1_2<8									
ormula input									
Variable		_	Operators	Functions					
BL1	*		+	and		*			
ita ta Plata OC									
ste to Flate QC									
Input data:	Diffe	erence d	lata			~			
Control 1	:	BL1	~	Mean:		s:			Ť
Control 2	2:	BL1	~	Mean:		s:			
Control 3	}:	BL1	~	Mean:		s:			
Control 4	k:	BL1	~	Mean:		s:			Ť
<ul> <li>All works</li> </ul>	spaces								
🔿 Workspa	aces create	d during	the last	days					
O Workspa	aces contai	ning nar	me						
	Validation C NC1_1<8 NC1_2<8 NC1_2<8 mula input Variable BL1 Late to Plate QC Input data: Control 1 Control 1 Control 2 Control 4 Control 4 Control 4	Validation Conditions NC1_1<8 NC1_2<8 Inc1_2<8 Input Variable BL1 Variable BL1 Control 1: Control 1: Control 2: Control 3: Control 3: Control 4: Other Spaces Workspaces create Vorkspaces create Vorkspaces contai	Validation Conditions         NC1_1<8         NC1_2<8         Input         Variable         BL1         Variable         BL1         Input data:         Difference d         Control 1:         BL1         Control 2:         BL1         Control 3:         BL1         Control 4:         BL1         Ontrol 3:         BL1         Outrol 4:         BL1	Validation Conditions         NC1_1<8         NC1_2<8         mula input         Variable       Operators         BL1       • • • •         te to Plate QC         Input data:       Difference data         Control 1:       BL1 •         Control 2:       BL1 •         Control 3:       BL1 •         Control 4:       BL1 •         Ocntrol 4:       BL1 •         Overkspaces       • All workspaces created during the last         Workspaces containing name       1	Validation Conditions         NC1_1<8         NC1_2<6         rmula input         Variable       Operators         BL1       Image: Control 1:         Control 1:       BL1         Control 2:       BL1         Control 3:       BL1         Control 4:       BL1         Variable       Mean:         Operators       Mean:         Control 1:       BL1         Mean:       Mean:         Operators       Image: Control 2:         BL1       Mean:         Operators       Image: Control 3:         BL1       Mean:         Operators       Image: Control 4:         Operators       Image: Control 4:         Norkspaces       Image: C	Validation Conditions         NC1_1<8         NC1_2<6         rmula input         Variable       Operators         BL1       Image: Control 1:         Input data:       Difference data         Control 1:       BL1         Control 2:       BL1         Control 3:       BL1         Control 4:       BL1         Ø       All workspaces         Workspaces containing name       1	Validation Conditions         NC1_1<8         NC1_2<8         rmula input         Variable       Operators         BL1       Image: Stress of the	Validation Conditions         NC1_1<8         NC1_2<8         rmula input         Variable       Operators         Functions         BL1       •         •       •         Input data:       Difference data         •       •         •       Mean:         •       •         •       Control 1:         BL1       Mean:         •       Control 3:         BL1       Mean:         •       Control 4:         BL1       Mean:         •       All workspaces         •       All workspaces containing name	Validation Conditions   NC1_1<8   NC1_2<8     Imula input     Variable   Operators   BL1   Imut data:     Difference data     Input data:   Difference data     Imut data:     Difference data     Imut data:   Difference data     Imput data:     Difference data     Imput data:   Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data:     Difference data     Imput data: <t< th=""></t<>



# 6.2.3 Organizarea raportului tipărit

În bara de control, dați clic pe **Printed repor**t (Raport tipărit) din item-ul **Data handling (Prelucrare date)**. Este afișat următorul ecran:

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       □       E       Data Collection       Up         Append       □       Image: Collection       Down         Vertical list       □       Image: Collection       Down         Image: Collection       □       Image: Collection       Image: Collection       Down         Image: Collection       □       Image: Collection       Image: Collection       Image: Collection       Down         Image: Collection       Image: Collection <t< th=""></t<>					
E Signature	Import					

Pe bara **Data selection (Selectare date)**, toate datele disponibile ale raportului se găsesc în caseta **Available data (Date disponibile)**. Utilizând butoanele **Insert and Append (Introducere şi Anexare)**, datele pot fi transferate în caseta **Selected data (Date selectate)**. Datele pot fi de asemenea transferate, utilizând funcția selectare şi introducere.

În caseta **Print as (Tipărire ca)**, alegeți între tipărirea datelor ca o matrice sau ca o listă cu o orientare specială.

În acest exemplu, trebuie creat un raport care să conțină parametrii de măsură, dispunere placă, valori blanked, curba standard, concentrațiile IgM, definiție valoare critică, rezultate calitative ale probelor și validărilor.

Înainte de crearea raportului, Vertical list/Difference data (Listă verticală/Date diferențiale) default trebuie îndepărtată din caseta Selected data (Date selectate). Numai Measurement parameters (Parametrii de măsură) mai rămân în caseta Selected data (Date selectate). Print as List (Tipărire ca listă) trebuie schimbat în Print as Matrix (Tipărire ca matrice).



Print as O 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:	
<ul> <li>Instrument data</li> <li>Reduced data</li> <li>Difference data</li> <li>Difference data - Mean</li> <li>Difference data - Mean</li> <li>Difference data - Variation co</li> <li>Transformed data</li> <li>Concentrations</li> <li>Qualitative Results</li> <li>'s Sample IDs</li> <li>Method Iayout</li> <li>Qualitative Results</li> <li>Remarks</li> <li>Error protocol</li> <li>Kendit Trail</li> <li>Signature</li> <li>Signature</li> </ul>	Insert  Append  Append	n es e all
	Import	

Selectați Method layout/Layout (Configurație metodă/Configurație) din caseta Available data (Date disponibile) și atașați-o la raport, dând clic pe Append (Atașament). Apoi introduceți Blank reduction (Reducere valoare blank), Mean conc. (UA/ml)(Conc. medie) și Cutoff results (Rezultate valori critice) în matrice selectând item-urile corespunzătoare și dând clic pe Insert (Introducere).

Append (Ataşament) Graph (Grafic): Standard curve (Curbă standard), Cutoff definition (Definiție valoare critică) și QC Validation criteria (Criterii Validare QC) la datele selectate. Partea de setare a datelor a procedurii de definire a raportului este completă; caseta de dialog Printed report (Raport tipărit) arată astfel:





### 6. Exemplu de aplicație

Pe barele **Header (Antet)** și **Footer (Subsol)**, definiți dispunerea antetului și subsolului raportului (a se vedea Manualul de utilizare Magellan pentru detalii suplimentare).

# Data Export (Export de date)

În bara de control, selectați **Data export (Export de date)** din item-ul **Data** handling (Prelucrare date). În acest exemplu, rezultatele de dispunere și separare trebuie stocate ca fișier ASCII. Selectați rezultatele Layout (Dispunere) și Cutoff (Valoare critică) din fereastra Available data (Date disponibile); dați clic pe  $\rightarrow$  săgeată pentru a le introduce în fereastra Selected data (Date selectate). Ecranul afișează următoarele informații:

Available data:			Selected data:	
⊕… 🛍 Instrument data	^		Layout	Up
🖨 🔆 Reduced data			Difference data	Dawn
Difference data				Down
Difference data - Mean				
Difference data - Standard deviation				
Difference data - Variation coefficie				
🕀 者 Transformed data				
🕀 頂 Concentrations		•		
Cutoff results				
⊕ <sup>1</sup> ², Sample IDs		<u></u>		
🖃 👜 Method layout				
Well positions				
Strip method names	-			
🔄 🔄 Original Concentrations				
Dilution factors				
	_			]
Export Options Export to ASCIL File	Г	Export t	o Excel	
	-	Liport		



Indicație Datele exportate trebuie să conțină întotdeauna Layout or Sample ID List (Lista ID-urilor de dispunere și de probe).



# Automated data handling (Prelucrarea automată a datelor)

În bara de control, selectați Automated data handling (Prelucrare automată a datelor) din item-ul Data handling (Prelucrare date).

Automatic	
📃 load sample ID list	More
save workspace	More
export to ASCII file	
export to ASTM file (LIS)	More
export to Excel	
🔲 print	
view results after measurement	More

Selectați export to ASCII file (export ca fișier ASCII), și view results after measurements (vizualizare rezultate după măsurări). În Magellan Tracker, save workspace (salvare spațiu de lucru) este selectat din setarea inițială și nu poate fi modificat.



# Salvarea metodei

Dați clic pe **Next (Continuare)** pentru a deschide fereastra **Save as (Salvare ca)**. Introduceți numele de fișier al metodei și completați orice alt câmp, dacă este necesar.

Create/Ed	lit a Method			
Save in:	🗁 mth	Name 🔺	Remarks	Status
Filenamer	Method1 mth			
File remarks	mediourtinut			
The remains	5.			
Audit trail c	omment:			
Org	anize favorites Audit trail	Signatures	Method password:	
L H	ncel	Ru	n this method now SAVE&FI	NISH
			© 2009	Tecan

Câmp text Filename (nume fişier)	Trebuie introdus un nume de fişier. Este sugerat un nume inițial de fişier în mod automat, dar acesta poate fi modificat.
Câmp text File remarks (Observaţii fişier)	Comentariile introduse aici vor fi salvate şi afişate cu numele fişierului.
Câmp text Comentariu Audit trail (Pistă de audit)	Comentariile introduse aici vor fi stocate în pista de audit. Această opțiune este disponibilă numai cu <b>Magellan Tracker</b> .
Butonul Organize Favorites… (Organizare favoriți)	Apare caseta de dialog <b>Organize Favorites</b> (Organizare favoriți).
Câmp text Method password (parolă metodă)	Introduceți o parolă a metodei pentru a proteja metoda.
Casetă de verificare Run this method now (Rulare acum a acestei metode)	Această metodă va rula imediat după ce se dă clic pe Save&Finish (Salvare și Terminare).

Pentru detalii suplimentare, vă rugăm să consultați Manualul de utilizare Magellan.



# 6.2.4 Run the Method (Rulare metodă)

Dacă este selectată Run this method now (Rulare acum a acestei metode) în caseta de dialog Save as (Salvare ca) a Create/edit a method wizard (Asistentului Creare/Editare metodă), caseta de dialog Start Measurement Wizard/Start Measurement (Asistent Start Măsurare/Start măsurare) va apărea după ce se dă clic pe Save (Salvare). În dialogul Start Measurement (Start măsurare) este alocat un nume inițial al spațiului de lucru și pote fi modificat de către utilizator, dacă se dorește acest lucru.

Start Measurement			
Measurement			Measurement parameters
Workspace: Method: Sample ID list: Instrument Use state Plate in Movements	05112009-001.wsp Method1.mth cker Temp. control	Arb. cycle kin Modify layout Insert Current: n. def °C Target: n. def °C Optimize Z-position	Plate Plate Description: [GRE96ft] - Greiner 96 Flat Tr- Plate with Cover: No Barcode: No Part of Plate Range: A1:H12 Absorbance Measurement wavelength: 492 nm Reference wavelength: 620 nm Label: Label1
Please note:	<<< Back		START
			© 2009 Tecan

Dați clic pe **Start (Start)** pentru a începe măsurarea. Va fi creat un spațiu de lucru în mod automat, care conține toate informațiile introduse anterior și va colecta toate valorile de măsură. În timp ce măsurare este executată, apare o casetă de dialog despre starea măsurării, indicând progresul măsurării.

După ce măsurarea s-a încheiat, apare caseta de dialog **Results (rezultate)**, în care pot fi vizualizate toate rezultatele și calculele. Mesajele de eroare pot apărea, atunci când se execută o măsurare fără lichidele corepunzătoare (de ex. standarde).



# 6.2.5 Evaluarea rezultatului

Selectați **Evaluate results (Evaluare rezultate)**pentru a vizualiza și a evalua datele brute. Parametrii de evaluare pot fi vizualizați și datele pot fi reevaluate.

Această secțiune vă ghidează prin **Evaluate Results wizard (Asistentul Evaluare rezultate)** utilizând un fișier exemplu de spațiu de lucru instalat după instalarea software-ului Magellan.

În caseta de dialog *Wizard List (Listă asistenți)*, dați clic pe **Evaluate results** (Evaluare rezultate).

Dați clic pe Next (Continuare) de pe pagina Welcome (de întâmpinare) a Evaluate Results wizard (asistentului Evaluare rezultate) și apare caseta de dialog Select a file (selectați un fișier).

Selectați spațiul de lucru Quantitative ELISA

**example\_Sunrise\_InfiniteF50.wsp** din lista fişierelor şi daţi clic pe **Make your selection (Selectaţi)**. Calculele sunt executate şi este afişată următoarea fereastră de dispunere a plăcii:

File Edit Instrument	el ISA e	xample_Sunn	se_Infiniter	50.wsp - Q	uantitative l	LISA examp	le_Sunrise.	nth							<u> </u>
EdR		10% 10	0% Zoor	n											
	\$														
Instrument data (>) Measurement data Reference data	Evaluate r	0	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	Edit method		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	
Reference data as Colors           Reduced data         Image: Colors           Difference data         Image: Colors	-	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 S		E	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	SM1_12 2/2	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	N Austria
Miscellaneous (A) Remarks Error protocol		H	ST1_3 1/2	SM1_1 1/2	SM1_5 1/2	SM1_9 1/2	SM1_13 1/2	SM1_17 1/2	SM1_21 1/2	SM1_25 1/2	SM1_29 1/2	SM1_33 1/2	SM1_37 1/2		TECA
Aut. rai Spritze															
	M ~ SI PI	leasurement para UNRISE late late Description: [	meters SUNRISE_Stan	dard] - Greiner	96 Flat										
Help Cancel <<< Back														NEXT	
													© 200	9 Tecan	

În fiecare godeu este afişată valoarea calculată. În funcție de item-ul selectat din bara de control, fereastra de dispunere a plăcii se modifică în mod corespunzător. Parametrii și setările pot fi modificate, utilizând item-urile din bara de control. Dacă metoda trebuie modificată, dați clic pe bara **Edit method** (Editare metodă).



Dați clic în godeu cu tasta dreaptă a mouse-ului și apare meniul sensibil la context:

SM1	_10 SM1_14 SM1_18
2 0.1	Summary Details Edit
SM <sup>2</sup> 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>2</sup>	Graph: Spectra Graph: Dilution series
0.0	Mask/Unmask selection Show/Hide layout

Selectând **Summary (Rezumat)** este afișată următoarea fereastră, asigurând informații detaliate despre definiția și setările godeului ales:

Well: C5					
🗩 📠 Method layout					
🕀 🛲 Identifier					
🕀 📠 Liquid Alias					
🗉 👜 Replicate					
🗉 📠 Dilution					
🗉 🛲 Multilabel data reduction					
🗉 🛲 Transformation formula(s)					
Concentration transformation formula(s)					
🕀 📠 Strip method names					
🕀 🛍 Instrument data	🛨 🛗 Instrument data				
🗄 💥 Reduced data					
🕀 🔟 Averages					
🕀 者 Transformed data					
🕀 🗊 Concentrations					
🗄 🔟 Statistics					
left Up right					
Expand All	ОК				

Dați clic pe **Next (Continuare)** în fereastra dispunerii plăcii și apare caseta de dialog **Save as (Salvare ca)**, unde puteți introduce un nume de fișier și observații. Dați clic pe butonul **Save (Salvare)** de pe partea stângă a ferestrei pentru a salva fișierul; puteți continua lucrul la metodă sau spațiul de lucru. Dați clic pe butonul **Finish (Terminare)** de pe partea dreaptă a ecranului pentru a salva fișierul și pentru a închide asistentul. Programul revine la lista asistenților.



# 6.2.6 Rezumatul definiției ELISA cantitative în Magellan

#### 1. Valoare blank de scădere

#### Definiții în Magellan

Dați clic pe Add new transformation (Adăugare transformare nouă) din bara de control și apare o fereastră, întrebându-vă dacă doriți să definiți o Blank reduction (Reducere valoare blank). Dați clic pe Yes (Da) și formula Blank reduction (Reducere valoare blank) este alocată în mod automat tuturor godeurilor.

#### 2. Definirea concentrațiilor

Definiții în Magellan (Control bar – Method Iayout/ Conc.-, Dil.-, Ref.values (Bară de control – Configurare metodă/ Valori conc., dil., ref.))

Identificator selectat: ST Unitate: UA/ml

ST1_1	5	(ST1_1Standard 1 primul grup experimental)
ST1_2	10	(ST1_2Standard 2 primul grup experimental)
ST1_3	20	(ST1_3Standard 3 primul grup experimental)
ST1_4	40	(ST1_4Standard 4 primul grup experimental)
ST1_5	80	(ST1_5Standard 5 primul grup experimental)
ST1_6 1	60	(ST1_6Standard 6 primul grup experimental)

## 3. Definirea curbei standard

# Definiții în Magellan (Control bar – Concentrations/ Standard curve (Bară de control - Concentrații/Curbă standard))

Input data (Date de intrare)	blank reduction (reducere valoare blank)
Analysis type (tip analiză)	linear regression (regresie liniară)
X-axis (ax X)	liniar
Y-axis (ax Y)	liniar

#### 4. Definire valori critice

# Definiții în Magellan (Control bar – Evaluate data/ Cutoff definition (Bară de control - Evaluare date/ Definire valoare critică))

Input data (Date de intrare):	Mean conc. (UA/ml) (Conc. medie)
Limite:	22
	18

Positive (Pozitiv)  $\ge$  22 > intermediate (intermediar)  $\ge$  18 > negative (negativ) Non-competitive test (Text necompetitiv)

#### 5. Validare QC

Definiții în Magellan (Control bar – Evaluate data/ QC validation (Bară de control - Evaluare date/ Validare QC)):

Input data (Date de intrare): Single conc. (UA/ml) (Conc. unică)

Condiție validare 1	NC1_1<8
---------------------	---------

Condiție validare 2 NC1\_2<8

NC1\_1.....Control negativ prima reproducere primul grup experimental

NC1\_2.....Control negativ a doua reproducere primul grup experimental



# 7. Curăţarea, întreţinerea şi eliminarea deşeurilor

# 7.1 Introducere

Acest capitol conține următoarele proceduri:

- Curățarea instrumentului
- Dezinfectarea instrumentului
- Întreținerea instrumentului
- Introducerea sau înlocuirea filtrelor în centrifuga cu filtre
- Instrucțiuni de eliminare a deșeurilor









# 

ÎNAINTEA CURĂŢĂRII ȘI DEZINFECTĂRII, DECONECTAȚI INSTRUMENTUL DE LA SURSA EXTERNĂ DE ENERGIE ELECTRICĂ.

#### PRECAUŢIE

ÎNDEPĂRTAȚI MANUAL DISPOZITIVUL DE TRANSPORT AL PLĂCII NUMAI DACĂ INSTRUMENTUL A FOST OPRIT.

# 7.2 Curăţarea instrumentului



# AVERTIZARE

PROCEDURA DE CURĂȚARE TREBUIE EXECUTATĂ ÎNTR-O CAMERĂ BINE AERISITĂ, DE CĂTRE PERSONAL AUTORIZAT CALIFICAT ECHIPAT CU MĂNUȘI RECICLABILE, PRECUM ȘI OCHELARI ȘI ÎMBRĂCĂMINTE DE PROTECȚIE.

Curăţaţi carcasa dispozitivului şi dispozitivul de transport al plăcii cu o cârpă uscată şi umezită. Dacă este foarte murdară, curăţaţi-o cu o cârpă umezită cu o soluţie de 70% etanol sau detergent delicat, Microcide SQ sau Decon 90. Ştergeţi până la uscare cu un prosop din hârtie.

Dacă instrumentul este stropit cu orice fel de lichid, acesta trebuie îndepărtat imediat pentru a preveni scurgerea lichidului în sistemul optic, cauzând pierderea performanței sau o eroare.



# 7.3 Dezinfectarea instrumentului



#### AVERTIZARE

#### DACĂ LICHIDUL VĂRSAT PE DISPOZITIVUL DE TRANSPORT AL PLĂCII ESTE POSIBIL INFECȚIOS, ACESTA VA FI DEZINFECTAT CONFORM LEGILOR ȘI REGLEMENTĂRILOR NAȚIONALE RELEVANTE.

Toate piesele instrumentului care intră în contact cu probe biologice, probe ale pacienților, probe pozitive de control sau material periculos, trebuie tratate ca zone posibil infecțioase.



# AVERTIZARE

PROCEDURA DE DEZINFECTARE ȘI DEZINFECTANȚII TREBUIE SĂ SE CONFORMEZE LEGILOR ȘI REGLEMENTĂRILOR NAȚIONALE RELEVANTE.



# AVERTIZARE

ESTE FOARTE IMPORTANT CA INSTRUMENTUL SĂ FIE DEZINFECTAT ÎNAINTE DE A FI MUTAT DIN LABORATOR SAU ÎNAINTEA EXECUTĂRII ORICĂREI LUCRĂRI CU ACESTA.

Înainte ca instrumentul să fie înapoiat reprezentantului local de vânzări sau unui centru de service, toate suprafeţele şi suportul plăcii trebuie dezinfectate şi trebuie completat un certificat de siguranţă de către autoritate operaţională. Dacă nu este asigurat un certificat de siguranţă, este posibil ca instrumentul să nu fie acceptat de către reprezentantul local de vânzări sau de centrul de service, sau autoritățile vamale îl pot reține.

# 7.3.1 Soluții de dezinfectare

Suprafețele exterioare și suportul plăcii instrumentului trebuie dezinfectate utilizând un dezinfectant, precum:

- Microcide SQ
- Decon 90
- 70 % etanol



AVERTIZARE PERICOL DE INCENDIU ȘI EXPLOZIE!

SOLUŢIILE DE ALCOOL, PRECUM ETANOL SAU IZOPROPAN, SUNT INFLAMABILE ȘI ATUNCI CÂND SUNT MANIPULATE ÎN MOD INADECVAT POT CAUZA EXPLOZII ȘI/SAU INCENDIU. TREBUIE RESPECTATE PRESCRIPȚIILE CORESPUNZĂTOARE DE LABORATOR PRIVIND SIGURANȚA.



PRECAUȚIE NU UTILIZAȚI ACETONĂ, DEOARECE DETERIOREAZĂ CAPACELE.



# 7.3.2 Procedura de dezinfectare

Dacă laboratorul nu deține nicio procedură specifică de dezinfectare, trebuie utilizată următoarea procedură pentru a dezinfecta suprafețele exterioare și suportul plăcii instrumentului.







# AVERTIZARE

PROCEDURA DE DEZINFECTARE TREBUIE EXECUTATĂ ÎNTR-O CAMERĂ BINE AERISITĂ, DE CĂTRE PERSONAL AUTORIZAT CALIFICAT ECHIPAT CU MĂNUŞI RECICLABILE, PRECUM ȘI OCHELARI ȘI ÎMBRĂCĂMINTE DE PROTECȚIE.

#### PRECAUŢIE

DEZINFECTANTUL DE SUPRAFEȚE POATE INFLUENȚA ÎN MOD NEGATIV PERFORMANȚA INSTRUMENTULUI DUMNEAVOASTRĂ, DACĂ ESTE APLICAT SAU DACĂ INTRĂ ÎN MOD ACCIDENTAL ÎN INSTRUMENT.

#### AVERTIZARE

ÎNAINTE DE DEZINFECTARE, DECONECTAȚI INSTRUMENTUL DE LA SURSA DE ALIMENTARE CU ENERGIE ELECTRICĂ, PENTRU A EVITA PERICOLUL UNUI INCENDIU SAU UNEI EXPLOZII.

- 1. Purtați mănuși, ochelari și îmbrăcăminte de protecție.
- 2. Pregătiți un coș adecvat pentru toate deșeurile rezultate în urma procedurii de dezinfectare.
- 3. Deconectați instrumentul de la software și opriți instrumentul.
- 4. Scoateți cu atenție suportul plăcii din instrument.



- Aplicaţi cu atenţie soluţia de dezinfectare pe suportul plăcii, conform Manualului de utilizare al producătorului. Nu utilizaţi prea mult dezinfectant, pentru a preveni scurgerea soluţiei în instrument sau murdărirea lentilelor atunci când suportul plăcii este introdus în instrument.
- După timpul necesar de contact (conform Manualului de utilizare al producătorului), ştergeţi suportul plăcii utilizând un prosop moale din hârtie umezit cu un detergent delicat sau apă distilată, pentru a îndepărta urmele de dezinfectant.



# 7. Curățarea, întreținerea și eliminarea deșeurilor

7. Introduceți cu atenție suportul plăcii în instrument.



- 8. Aplicați cu atenție dezinfectantul pe placa de bază a suportului plăcii.
- 9. După timpul necesar de contact, ștergeți placa de bază a suportului plăcii, utilizând un prosop moale din hârtie umezit cu un detergent delicat sau apă distilată, pentru a îndepărta urmele de dezinfectant
- 10. Aplicați cu atenție soluția de dezinfectare pe toate suprafețele exterioare ale instrumentului.
- 11.După timpul necesar de contact, ştergeţi instrumentul, utilizând un prosop moale din hârtie umezit cu un detergent delicat sau apă distilată, pentru a îndepărta toate urmele de dezinfectant.
- 12.Uscați prin ștergere suprafața instrumentului cu un prosop moale din hârtie.
- 13.Repetați procedura de dezinfectare pe orice accesorii, care sunt mutate sau înapoiate.
- 14. Aruncați coșul cu deșeuri, conform legilor și reglementărilor naționale relevante.
- 15.Dezinfectați-vă mâinile și curățați-le cu un detergent delicat.

Atunci când înapoiați instrumentul reprezentantului local de vânzări/centrului de service, continuați cu următoarele etape:

16.Împachetați instrumentul și accesoriile sale.

17.Completați certificatul de siguranță (a se vedea mai jos) și atașați-l la partea exterioară a cutiei, astfel încât să fie vizibil explicit.

# 7.3.3 Certificat de siguranță

Pentru a asigura siguranța și sănătatea personalului, clienții noștri sunt rugați să completeze un **Certificat de siguranță** (care a fost furnizat cu instrumentul) și să atașeze o copie la partea superioară a containerului în care este returnat instrumentul (vizibil de la partea superioară a containerului de transport!) și o altă copie la documentele de transport, înainte de a-l expedia centrului de service sau reparație.

Instrumentul trebuie dezinfectat la sediul autorității operaționale înainte de expediere (a se vedea 7.3.2 Procedura de dezinfectare).

Procedura de dezinfectare trebuie executată într-o cameră bine aerisită, de către personal autorizat și calificat, echipat cu mănuși reciclabile, fără pudră, ochelari și îmbrăcăminte de protecție.

Procedura de dezinfectare trebuie executată conform reglementărilor naţionale, regionale și locale.



Dacă nu este asigurat un Certificat de siguranță, este posibil ca instrumentul să nu fie acceptat de centrul de service.

Serviciul dvs. local de asistență clienți Tecan vă poate trimite o nouă copie a Certificatului de siguranță, dacă este necesar.

# 7.4 Plan de întreţinere preventivă pentru INFINITE F50 PLUS

Următoarele proceduri de întreținere preventivă sunt recomandate.

# 7.4.1 Lunar

Curățați carcasa și dispozitivul de transport al plăcii cu un detergent delicat cel puțin o dată pe an; mai des, dacă este necesar.



PRECAUȚIE NU UTILIZAȚI ACETONĂ, DEOARECE DETERIOREAZĂ CAPACELE.

# 7.4.2 La fiecare 4 ani

Se recomandă înlocuirea filtrelor la fiecare 4 ani.

# 7.5 Înlocuirea filtrelor și instalarea

INFINITE F50 PLUS trebuie conectat la Magellan, pentru a realiza o introducere sau o înlocuire a filtrelor, ghidată de software. Dacă conexiunea se întrerupe în timpul procedurii, din cauza întreruperii accidentale dintre instrument și calculator, Magellan trebuie închis și instrumentul trebuie oprit. În acest caz, continuați procedura, așa cum se indică mai jos. După terminare, restabiliți conexiunea prin repornirea dispozitivului și a software-ului Magellan, și definiți filtrele nou introduse.



# PRECAUŢIE

ATUNCI CÂND MANIPULAȚI FILTRELE, FIȚI ATENȚI SĂ NU LE ZGÂRIAȚI SAU SĂ NU LE MURDĂRIȚI CU DEGETELE SAU CU PRAF.



# 7.5.1 Procedura de schimbare a filtrelor

Filtrele centrifugei standard cu filtre pot fi înlocuite sau suplimentate, utilizând următoarea procedură:

- 1. În fereastra Wizard list (Listă asistenți), dați clic pe Miscellaneous (Diverse).
- 2. Dați clic pe Instrument control (Control instrument)
- 3. Dați clic pe Define filter slides... (Definire ghidaje filtre)
- 4. Dați clic pe Filterswitching (Schimbare filtre) pentru a începe procedura.
- 5. Îndepărtați orice microplacă din suportul plăcii!
- 6. Înclinați instrumentul cu atenție spre spate, până când stă aşezat pe spate cu partea inferioară îndreptată spre dumneavoastră.
- Îndepărtaţi placa de acoperire de pe fundul instrumentului îndepărtând şuruburile portocalii.



8. Îndepărtați centrifuga cu filtre fixată magnetic, scoțându-o din instrument.



9. Aşezaţi centrifuga cu filtre pe o suprafaţă curată şi netedă.



# 7. Curățarea, întreținerea și eliminarea deșeurilor

10.Atunci când înlocuiți un filtru, utilizați unealta de asamblare a filtrului pentru a îndepărta filtrul din canalul pentru filtru.

Vă rugăm să contactați serviciul dvs. local de asistență clienți Tecan pentru unealta de asamblare a filtrului și pentru filtre disponibile.



- 11.Aliniați unealta de asamblare a filtrului cu crestătura inelului de oprire. Răsuciți unealta și îndepărtați inelul de oprire, trăgându-l afară din canalul pentru filtru.
- 12.Răsuciți centrifuga cu filtre, astfel încât filtrul să alunece în afară din canal. Nu utilizați unealta de asamblare a filtrului pentru a împinge filtrele în afară din canalul pentru filtru, deoarece filtrul poate fi zgâriat.
- 13. Trebuie introdus un nou filtru în canalul pentru filtru, în direcția corectă, având grijă să nu zgâriați filtrul sau să lăsați amprente pe acesta.





Indicație Asigurați-vă că filtrul este introdus corect.



# 7. Curăţarea, întreținerea și eliminarea deșeurilor

14.Puneți inelul de oprire pe capătul uneltei de asamblare a filtrului și răsuciți-l, astfel încât să nu poată aluneca în afară.



- 15.Utilizând unealta de asamblare a filtrului, împingeți inelul de oprire în canalul pentru filtru și apăsați-l la loc în mod ferm.
- 16.Răsuciți unealta, până când crestătura din inelul de oprire este aliniată cu capătul uneltei de asamblare a filtrului și îndepărtați unealta.



- 17.Puneți la loc centrifuga cu filtre în canalul centrifugei cu filtre și împingeți-o adânc până când este asigurată magnetic.
- 18.Remontați placa de acoperire pe fundul instrumentului cu cele trei şuruburi portocalii.
- 19. Aduceți la loc dispozitivul în poziție orientat în sus.
- 20.Dați clic pe **OK** pentru a încheia procedura și pentru a inițializa centrifuga cu filtre.
- 21.Definiți filtru nou introdus (a se vedea capitolul următor pentru procedura detaliată).


### 7.5.2 Definirea filtrelor

În caseta de dialog **Filter Definițion (Definiție filtru)** alocați lungimile adecvate de unde filtrelor înlocuite, introducând noile lungimi de undă în pozițiile corespunzătoare.

Dacă un filtru a fost introdus într-o poziție de filtru nou, activați poziția corespunzătoare a filtrului, selectând caseta de verificare și introducând lungime adecvată de undă.

Dând clic pe **Save (Salvare)**, definițiile de filtru vor fi salvate și filtrele vor fi inițializate.

De îndată ce filtrele au fost inițializate, instrumentul este pregătit pentru măsurări.



Indicație Aveți grijă să nu amestecați pozițiile filtrelor și lungimile de undă ale filtrelor, deoarece acest lucru va cauza date greșite de măsură.

### 7.6 Eliminarea deşeurilor

### 7.6.1 Introducere

Respectați procedurile de laborator pentru eliminarea deșeurilor biologice periculoase conform reglementărilor naționale și locale.

Acest capitol asigură instrucțiuni despre eliminarea legală a deşeurilor acumulate în urma utilizării INFINITE F50 PLUS.



### PRECAUȚIE RESPECTAȚI TOATE REGLEMENTĂRILE FEDERALE, NAȚIONALE ȘI LOCALE DE MEDIU ÎNCONJURĂTOR.

### 7.6.2 Eliminarea ambalajului

Ambalajul este fabricat din material reciclabil. Dacă nu intenționați să îl păstrați pentru utilizarea viitoare, de ex. pentru transport și depozitare, vă rugăm să eliminați ambalajul conform reglementărilor locale.

### 7.6.3 Eliminarea materialelor de lucru

### AVERTIZARE

PERICOLELE BIOLOGICE POT FI ASOCIATE CU DEȘEURILE (MICROPLACĂ) RULĂRII PROCESULUI PE INFINITE F50 PLUS ABSORBANCE READER.



TRATAȚI MIROPLACA UTILIZATĂ, ALTE DEȘEURI ȘI TOATE SUBSTANȚELE UTILIZATE ÎN CONFORMITATE CU DIRECTIVELE DE BUNĂ PRACTICĂ DE LABORATOR.

INTERESAȚI-VĂ DE PUNCTELE CORESPUNZĂTOARE DE COLECTARE ȘI METODELE APROBATE DE ELIMINARE A DEȘEURILOR DIN ȚARA SAU REGIUNEA DUMNEAVOASTRĂ.



### 7. Curățarea, întreținerea și eliminarea deșeurilor

### 7.6.4 Eliminarea instrumentului

Dacă aveți întrebări referitor la eliminarea dispozitivului, vă rugăm să contactați serviciul dvs. local de asistență clienți Tecan.

Grad de poluare	2 (IEC/EN 61010-1)
Metodă de eliminare	Deşeuri contaminate





DEZAFECTATE (WEEE) INFLUENŢELE NEGATIVE ASUPRA MEDIULUI ÎNCONJURĂTOR ASOCIATE CU TRATAREA DEŞEURILOR.

**ATENŢIE** DIRECTIVA 2012/19/UE DESPRE ECHIPAMENTELE ELECTRICE ȘI ELECTRONICE

- NU ARUNCAȚI ECHIPAMENTUL ELECTRIC ȘI ELECTRONIC CA DEȘEURI MENAJERE NESORTATE.
- COLECTAȚI SEPARAT ECHIPAMENTELE ELECTRICE ȘI ELECTRONICE SCOASE DIN FUNCȚIONARE.

### AVERTIZARE

ÎN FUNCȚIE DE APLICAȚII, ESTE POSIBIL CA PIESELE DISPOZITIVULUI INFINITE F50 PLUS SĂ FI INTRAT ÎN CONTACT CU MATERIAL BIOLOGIC PERICULOS.

- ASIGURAŢI-VĂ CĂ TRATAŢI ACEST MATERIAL CONFORM STANDARDELOR ŞI REGLEMENTĂRILOR ÎN VIGOARE PRIVIND SIGURANŢA.
- DECONTAMNAȚI TOATE PIESELE ÎNAINTE DE ELIMINARE.

Pentru informații suplimentare despre produs, vă rugăm să contactați:

Tecan Austria GmbH

Untersbergstrasse 1A

A-5082 Grödig/Salzburg

AUSTRIA/EUROPA

T +43 6246 8933 444

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

www.tecan.com

# 8. Remediere defecțiuni

### 8.1 Introducere

Microprocesorul central controlează și verifică funcțiile electronice, precum și măsurările, operațiile și rezultatele. Dacă microprocesorul detectează o eroare sau o procedură incorectă de operare, este afișat un mesaj de eroare pe calculator.

### 8.1.1 Tabel al mesajelor de eroare și al remedierii defecțiunilor

Următorul tabel oferă o scurtă descriere a mesajelor de eroare și acțiunile de remediere a defecțiunilor.



Indicație Dacă apar alte mesaje de eroare care nu sunt menționate în tabelul de mai jos, vă rugăm să contactați reprezentantul dumneavoastră local al serviciului de asistență clienți Tecan.

Mesaj de eroare	Descriere	Remediere defecțiuni
Eroare de sistem		
"Lid Open Error" (Eroare capac deschis)	Capac deschis la începerea măsurării	Închideți capacul și începeți din nou măsurare
"MTP Init Error" (Eroare inițializare MTP)	Suportul MTP nu a putut fi iniţializat	Problemă harware: Defecțiune electronică, curea ruptă sau suport MTP blocat mecanic
"MTP lost steps abs( <i>steploss</i> ) > <i>max_steploss</i> "	Etape MTP pierdute în timpul măsurării Steploss: numărul etapelor pierdute max_steploss: numărul etapelor pierdute permise	Problemă harware: Defecțiune electronică, mişcare grea din punct de vedere mecanic
"Filter lost steps abs( <i>steploss</i> ) > <i>max_steploss</i> "	Etape pierdute ale centrifugei cu filtre în timpul măsurării. Steploss: numărul etapelor pierdute max_steploss: numărul etapelor pierdute permise	Problemă harware: Piesă electronică defectă, mişcare grea din punct de vedere mecanic
"USB timeout"	Timeout în comunicația USB	Eroare de sistem – raportați la serviciul de asistență clienți
"Lamp Low! Minim: <i>minimum</i> , Maxim: <i>maximum</i> " Acest mesaj apare la versiunile de firmware până la V1.11.	Intensitatea măsurată a luminii nu a atins nivelul aşteptat dintre <i>minimum (minim)</i> și <i>maximum</i> <i>(maxim)</i>	Problemă harware: Defecțiune electronică, fibră ruptă
Începând cu versiunile de firmware V1.12 și superioare, mesajul "Lamp Low" este înlocuit cu: "Prepare REF check"	Intensitatea măsurată a luminii nu a atins nivelul așteptat dintre <i>minimum (minim)</i> și <i>maximum</i> <i>(maxim).</i>	Problemă hardware: defecțiune electronică, fibră ruptă



### 8. Remediere defecțiuni

Mesaj de eroare	Descriere	Remediere defecțiuni
Eroare de sistem		
"Wavelength Not Available ! (Lungime de undă indisponibilă!) Lungime de undă: <i>wavelength</i> nm"	Filtrul cu lungimea de undă <i>wavelength</i> nu a putut fi găsit în centrifuga cu filtre	Eroare de sistem – raportaţi la serviciul de asistenţă clienţi
"Channel Low! Canal: <i>channel_nr</i> , Minim: <i>minimum</i> , Maxim <i>maximum</i>	Semnalul de pe canalul <i>channel_nr</i> nu a atins zona aşteptată dintre <i>minimum</i> ( <i>minim</i> ) și <i>maximum (maxim</i> )	Problemă harware: Defecțiune electronică, fibră ruptă
"Invalid Wavelength! (Lungime de undă invalidă) Lungime de undă: <i>wavelength</i> nm"	Lungimea de undă a filtrului se află în afara domeniului de lungimi de undă ale LED-ului alb și albastru	Eroare de sistem – raportați la serviciul de asistență clienți
"Lamp Overflow!" Minim: <i>minimum</i> , Maxim: <i>maximum.</i>	Semnalul de pe ADC depăşeşte zona aşteptată dintre <i>minimum</i> <i>(minim)</i> şi <i>maxim</i>	Problemă harware: Defecțiune electronică
"Value Not Set: (Valoare nesetată) <i>value -1</i> "	Valoarea <i>value</i> nu este setată	Eroare de sistem – raportați la serviciul de asistență clienți
"Filter Init Error" (Eroare inițializare filtru)	Suportul filtrelor nu a putut fi iniţializat	Problemă harware: Defecțiune electronică, suport al centrifugei cu filtre blocat mecanic

### 8.1.2 Definiția 'Overflow'

Dacă rezultatul măsurării absorbanţei nu se află în cadrul specificaţiilor instrumentului (> 4.0 OD), va apărea un preaplin şi valoarea OD măsurată a godeului respectiv va fi înlocuită de 'Overflow'. Acest lucru se realizează de către software-ul de control şi nu de instrumentul în sine.

### 8.1.3 Cădere a tensiunii

În cazul căderii tensiunii, are loc următorul lucru:

Căderea tensiunii instrumentului, dar nu a calculatorului de control (de ex. calculator conectat la sistem de alimentare cu energie electrică, care nu poate fi întreruptă): Conexiunea USB între instrument și calculator se va întrerupe. Software-ul de control creează un mesaj de eroare.

Căderea tensiunii atât a instrumentului, cât și a calculatorului de control: calculatorul trebuie să se rebooteze. Nu va fi disponibilă nicio dată de măsură.

# 9. Abrevieri, mărci înregistrate şi simboluri

## 9.1 Abrevieri

Următoarele abrevieri sunt asigurate ca referință și pot apărea în Manualul de utilizare.

A	Amper
AC	Curent alternativ
ADC	Convertizor analogic digital
ANSI/SBS	Institutul American al Standardelor Naționale/Societatea pentru Ecranare Biomoleculară
ASCII	Codul Standardelor Americane pentru Schimb de Informații
ASTM	Societatea Americană pentru Testare și Material
°C	Grad Celsius
CE	Marcaj CE de conformitate
CFR	Codul Reglementărilor Federale
cm	Centimetru
DC	Curent direct
EC	Comunitatea Europeană
ELISA	Analiza pe bază de imunosorbenți enzimatici
EN	Norma Europeană
°F	Grade Fahrenheit
FDA	Administrarea alimentelor și medicamentelor
Hz	Hertz
IEC	Comisia Internațională Electrotehnică
ID	Identificare
IFU	Manual de utilizare
IQ	Calificarea instalării
IVD	Diagnostic in vitro
IVDR	Regulamentul privind diagnosticul in vitro (IVDR) (UE) 2017/746
kg	Kilogram
I	Litru
LED	Diodă electroluminescentă
LIS	Sistem Informatic de Laborator
mg	Miligram
ml	Mililitru
mm	Milimetru
MTP	Microplacă



### 9. Abrevieri, mărci înregistrate și simboluri

μΙ	Microlitru
NFM	Metal neferos
NIST	Institutul Național al Standardelor și Tehnologiei
nm	Nanometru
NRTL	Laborator de Testare Recunoscut Național
OD	Densitate optică
OQ	Calificare operațională
PCB	Printed Circuit Board (plăcii cu circuite imprimate)
RF	Frecvență radio
RoHS	Restricție pentru Utilizarea unor Anumite Substanțe Periculoase
SOP	Procedură Standard de Operare
USB	Bus Serial Universal
UA	Unități arbitrare
TÜV	Technischer Überwachungsverein (Asociația de Inspecție Tehnică)
V	Volt
VA	Volt Amper
WEEE	Echipamente electrice și electronice dezafectate

### 9.2 Mărci înregistrate

Următoarele nume de produse și orice mărci înregistrate și neînregistrate menționate în acest document sunt utilizate numai în scopuri de identificare și rămân proprietatea exclusivă a proprietarilor respectivi:

- Magellan<sup>™</sup>, Infinite<sup>®</sup>, MultiCheck<sup>™</sup>, Tecan<sup>®</sup> şi logoul Tecan Logo sunt mărci înregistrate ale Tecan Group Ltd., Männedorf, Elveţia
- Windows® şi Excel® sunt mărci înregistrate ale Microsoft Corporation, Redmond, WA, SUA
- Pentium® şi Atom<sup>™</sup> sunt mărci înregistrate ale Intel Corporation, Santa Clara, CA, SUA
- Adobe® Reader® este o marcă înregistrată a Adobe Systems Incorporated, Seattle, WA, SUA
- Microcide SQ<sup>TM</sup> este o marcă înregistrată a Global Biotechnologies Inc., Portland, ME, SUA
- Decon 90<sup>™</sup> este o marcă înregistrată a Decon Laboratories Ltd., Hove, East Sussex, UK



### 9.3 Simboluri

Următoarele simboluri apar pe instrument.

	Producător
$\sim$	Data fabricației
CE	Marcajul de conformitate CE
UK CA	Marcajul de evaluare a conformității pentru Regatul Unit arată că produsul etichetat respectă reglementarea aplicabilă în Marea Britanie.
i	Citiți Manualul de utilizare înainte de utilizarea instrumentului
IVD	Dispozitiv medical pentru diagnostic in vitro
UDI	Identificator unic al dispozitivului Simbolul UDI identifică suportul de date de pe etichetă.
REF	Număr comandă
SN	Numărul de serie
€ - - - - - - - - - - - - -	Etichetă USB
	Simbol WEEE
	Simbol RoHS, China
	NRTL TÜV SÜD MARK
	Periculos din punct de vedere biologic



# Index

### Α

Abrevieri	77
Absorbance (Absorbanță)	23
Asistent	
Ataşare semnătură	20
Creare/Editare metodă	20
Evaluare rezultate	19
Listă asistenți	19
Start măsurare	19
Asistentul	
Creare/Editare listă ID de probe	20

### В

### Bară de control

Acţiuni	
Cinetic	
Diverse	25
Măsurări	
Material de laborator	22
Bară de control Acțiuni	23
Bară de control Cinetic	24
Bară de control Diverse	25
Bară de control Măsurări	23
Bară de control Material de laborator	22

### С

Căderea tensiunii	76
Calificare operațională	39
Test MultiCheck	39
Testarea linearității	40
Testarea microplăcii	39
Testarea preciziei	40
Centrifugă cu filtre	35
Cerințe de energie	14
Cerințe de mediu înconjurător	14
Cerințe de sistem	15
Certificat de siguranță	68
Comment (Comentariu)	24
Control al calității	39
Control Bar (Bară de control)	22
Curățare	65

### D

Definirea filtrelor	73
Definirea măsurărilor	
Despachetarea	
Despachetarea și verificarea	
Procedura de despachetare	
Dezinfectare	65, 66
Certificat de siguranță	
Procedura	67
Soluții	
_	

### Ε

Editor parametrii de	măsură	21
----------------------	--------	----

Elemente de program	
(Timer) Aşteptare	25
Absorbanță	23
Agitare	23
Ciclu cinetic	24
Comentariu	24
Condiție cinetică	24
Incubare	
Indentare și Declanșare	
Parte a plăcii	
Placă	
Solicitare utilizator	
Eliminare	
Ambalaj	73
Instrument	74
Material de lucru	73
Eliminarea deseurilor	65, 73
Exemplu de aplicatie	,
Analiză cantitativă ELISA	
Exemplu de aplicatie	43

### I

Incubation (Incubare)	25
Info Pane (Panou Info)	27
Înlocuirea filtrelor	69
Instrument	
Accesorii	
Amplasare	
Caracteristici	
Descriere	
Specificații	
Întreținere	65, 69
Plan de întreținere	69

### Κ

Kinetic Condition (Condiție cinetică)	. 24
Kinetic Cycle (Ciclu cinetic)	. 24

### L

Listă	asistenți	19
-------	-----------	----

### М

Magellan	17
Magellan Tracker	17
Măgurări ciratica	20
	29
Másurári endpoint	27
Măsurări multilabel	28
Mesaje de eroare	75
Microplăci	37
Optimizarea performanţei	. 32

### 0

Optimizarea performanței 32	)
-----------------------------	---

### Ρ

Part of Plate (Parte a plăcii) ...... 22



### Index

Plate (Placă)	
PORNIREA instrumentului	
Preaplin	76
Procedura de auto-verificare	
Profilul utilizatorului	11
_	

### R

Remediere defecțiuni	75
-	

### S

Schimbarea filtrelor	70
Shaking (Agitare)	
Siguranța	7
Simboluri	79

Software 1	7
Calificarea instalării1	18
Instalarea 1	17
Interfaţa utilizatorului1	19
Listă asistenți 1	19
Specificații 3	36
U	
User Request (Solicitare utilizator)2	25
W	
Wait (Timer) ((Timer) Aşteptare) 2	25
Workflow Pane (Panou workflow) 2	26



# Serviciul de asistență clienți Tecan

Pentru întrebări sau asistență tehnică privind produsul dvs. Tecan contactați Serviciul local de asistență clienți Tecan. Pentru informații de contact accesați <u>http://www.tecan.com/</u>.

Înaintea contactării firmei Tecan pentru asistență privind produsul pregătiți următoarele informații în vederea asigurării unei asistențe tehnice optime (a se vedea plăcuța cu denumirea):

- Denumirea modelului produsului dvs.
- Numărul de serie (SN) al produsului dvs.
- Software și versiunea de software (dacă este cazul)
- Descrierea problemei și persoana de contact
- Data și ora apariției problemei
- Etapele deja parcurse pentru remedierea problemei
- Informațiile dvs. de contact (număr de telefon, număr de fax, adresă de e-mail etc.)

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:	Microplate Reader
Model Designation:	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.