# **•**TECAN•

Notice d'utilisation du lecteur d'absorbance

# INFINITE F50 PLUS et du logiciel MAGELLAN



Document n° :	30186914
	2023-06
N° de révision de document :	1.2
Version du micrologiciel :	3.33
Version du logiciel Magellan :	7.5







#### AVERTISSEMENT

#### LIRE ET SUIVRE ATTENTIVEMENT LES INSTRUCTIONS FOURNIES DANS CE MANUEL AVANT DE FAIRE FONCTIONNER L'APPAREIL.

#### Remarque

Tous les efforts ont été fournis afin d'éviter que des erreurs ne figurent dans le texte et les schémas ; cependant, Tecan Austria GmbH décline toute responsabilité en cas d'erreurs éventuelles présentes dans ce manuel.

La politique de Tecan Austria GmbH est de perfectionner les produits à mesure que de nouvelles techniques et de nouveaux composants arrivent sur le marché. Tecan Austria GmbH se réserve donc le droit de modifier les spécifications à tout moment *après vérification, validation, et obtention des approbations appropriées.* 

Tout commentaire sur ce manuel sera le bienvenu.



#### Fabricant

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

#### Copyright

Le contenu de ce manuel est la propriété de Tecan Austria GmbH. Toute copie, reproduction ou cession à des tiers est interdite sans autorisation écrite préalable.

Copyright © Tecan Austria GmbH Tous droits réservés. Imprimé en Autriche.

#### Déclaration pour certification UE

Voir la dernière page de la présente notice d'utilisation.

#### À propos de ce manuel

Notice originale. Ce document est prévu comme **Notice d'utilisation** pour le lecteur d'absorbance INFINITE F50 PLUS, qui est conçu pour mesurer l'absorbance lumineuse (densité optique) d'échantillons dans des microplaques à 96 puits. Elle doit servir de référence et de notice à l'utilisateur.

Ce document explique comment :

- Installer l'instrument
- Utiliser l'instrument
- Nettoyer et assurer la maintenance de l'instrument

#### Remarque sur les captures d'écran

Le numéro de version affiché dans les captures d'écran peut ne pas correspondre à celui de la version actuellement publiée. Les captures d'écran n'ont été remplacées que si le contenu relatif à l'application a changé.

#### Historique des modifications de la Notice d'utilisation

Version	Modifications
V 1.2	Nouvelle DoC (Declaration of Conformity)



# Avertissements, précautions et notes

Les types d'avertissements suivants sont utilisés dans ce manuel pour mettre en évidence des informations importantes ou pour avertir l'utilisateur de situations susceptibles d'être dangereuses :













PRÉCAUTION INDIQUE LA POSSIBILITÉ D'UNE DÉTÉRIORATION DU MATÉRIEL OU D'UNE PERTE DE DONNÉES SI LES INSTRUCTIONS NE SONT PAS SUIVIES.

#### **AVERTISSEMENT**

INDIQUE LA POSSIBILITÉ DE DOMMAGES CORPORELS, DANGER DE MORT OU DÉTÉRIORATION DU MATÉRIEL SI LES INSTRUCTIONS NE SONT PAS SUIVIES.

#### **AVERTISSEMENT**

INDIQUE LA PRÉSENCE POSSIBLE DE MATÉRIAUX PRÉSENTANT UN RISQUE BIOLOGIQUE. DES PRÉCAUTIONS DE SÉCURITÉ DE LABORATOIRE ADAPTÉES DOIVENT ÊTRE SUIVIES.

#### **AVERTISSEMENT**

CE SYMBOLE INDIQUE LA PRÉSENCE DE MATÉRIAUX INFLAMMABLES OU D'UN RISQUE D'INCENDIE. DES PRÉCAUTIONS DE SÉCURITÉ DE LABORATOIRE ADAPTÉES DOIVENT ÊTRE SUIVIES.

#### ATTENTION

DIRECTIVE 2012/19/UE SUR LES DÉCHETS D'ÉQUIPEMENTS ÉLECTRIQUES ET ÉLECTRONIQUES (DEEE)

EFFETS NÉGATIFS SUR L'ENVIRONNEMENT ASSOCIES AU TRAITEMENT DES DÉCHETS ;

- NE TRAITEZ PAS LES DÉCHETS D'ÉQUIPEMENTS ÉLECTRIQUES ET ÉLECTRONIQUES COMME DES DÉCHETS MÉNAGERS NON-TRIES.
- COLLECTEZ SÉPARÉMENT LES DÉCHETS D'ÉQUIPEMENTS ÉLECTRIQUES ET ÉLECTRONIQUES.



# Table des matières

1.	Séc	urité		7
	1.1	Sécurit	té de l'instrument	7
2.	Gén	éral		9
	2.1	Utilisat	tion prévue	
	2.2	Princip	pe de fonctionnement	
	2.3	Profils	d'utilisateur	
		2.3.1	Utilisateur professionnel – Niveau Administrateur	
		2.3.2	Utilisateur final ou utilisateur routinier	
		2.3.3	Techniciens de maintenance	
3.	Pris	e en mai	in	13
	3.1	Déballa	age et inspection	13
	•	3.1.1	Inspection de l'emballage livré	13
		3.1.2	Procédure de déballage	
	3.2	Caract	éristiques d'alimentation nécessaires	
	3.3	Caract	éristiques environnementales nécessaires	14
	3.4	Caract	éristiques système nécessaires	15
	3.5	Mettre	l'instrument sous tension	
	3.6	Logicie	el	17
		3.6.1	Introduction/ Utilisation prévue	17
		3.6.2	Procédure d'installation	17
		3.6.3	Qualification de l'installation	
		3.6.4	Commencer à travailler avec Magellan	
	3.7	Magell	an - Éditeur de paramètres de mesure	21
		3.7.1	Barre de contrôle	22
		3.7.2	Volet Flux de travail	
		3.7.3	Volet d'info	
	3.8	Magell	an - Définir les mesures	27
		3.8.1	Définir les mesures de point final	27
		3.8.2	Définir des mesures multi-étiquetage	
		3.8.3	Définir des mesures cinétiques	29
		3.8.4	Indentation et libération des éléments de programmes	
	3.9	Optimi	sation des performances	
		3.9.1	Emplacement de l'instrument	
		3.9.2	Procédure opérationnelle	
4.	Fon	ctionnal	ités de l'instrument	33
	4.1	Fonctio	onnalités de l'instrument	33
		4.1.1	Agitation de la microplaque	
	4.2	Descri	ption de l'instrument	33
	4.3	Descri	ption de la roue porte-filtres	35
	4.4	Spécifi	ications de l'instrument	
		4.4.1	Spécifications générales	
		4.4.2	Spécifications des mesures	
		4.4.3	Microplaques	
	4.5	Access	soires de l'instrument	
5.	Con	trôle qu	alité	39
	5.1	Introdu	uction	39
	5.2	Procéd	dure d'auto-vérification	39
	5.3	Qualifi	cation opérationnelle (QO)	39
		5.3.1	Test MultiCheck	
		5.3.2	Test de microplaque	
		5.3.3	Précision de l'instrument avec des échantillons liquides	40
		5.3.4	Linéarité de l'instrument avec des échantillons liquides	41
6.	Exe	mple d'a	application	43
	6.1	Introdu	uction	43
	6.2	Exemp	ble étape par étape : ELISA quantitatif	43



6.2.2       Créer une méthode       4         6.2.3       Organiser le rapport imprimé       5         6.2.4       Exécuter la méthode       5         6.2.5       Évaluer le résultat       6         6.2.6       Résumé de la définition de l'ELISA quantitatif dans Magellan       6
6.2.3       Organiser le rapport imprimé
6.2.4       Exécuter la méthode       50         6.2.5       Évaluer le résultat       60         6.2.6       Résumé de la définition de l'ELISA quantitatif dans Magellan       60
6.2.5       Évaluer le résultat
6.2.6 Résumé de la définition de l'ELISA quantitatif dans Magellan62
7. Nettoyage, maintenance et élimination des déchets6
7.1 Introduction
7.2 Nettoyer l'instrument
7.3 Désinfection de l'instrument
7.3.1 Solutions de désinfection
7.3.2 Procédure de désinfection
7.3.3 Certificat de sécurité
7.4 Plan de maintenance préventive pour INFINITE F50 PLUS6
7.4.1 Tous les mois
7.4.2 Tous les quatre ans
7.5 Remplacement et installation du filtre 6
7.5.1 Procédure de changement de filtre66
7.5.2 Définir les filtres7
7.6 Élimination des déchets
7.6.1 Introduction
7.6.2 Élimination du matériel d'emballage7
7.6.3 Élimination du matériel de fonctionnement7
7.6.4 Élimination de l'instrument
8. Dépannage
8.1 Introduction
8.1.1 Tableau des messages d'erreur et de dépannage
8.1.2 Définition de 'Overflow' (débordement)
8.1.3 Panne électrique
9. Abréviations, margues et symboles
9.1 Abréviations
9.2 Margues de commerce
9.3 Symboles
Index
Service client Tecan



# 1. Sécurité

# 1.1 Sécurité de l'instrument

- 1. Respectez toujours les précautions de sécurité de base lors de l'utilisation de ce produit, afin de réduire les risques de blessure, d'incendie ou de choc électrique.
- 2. Vous devez lire et comprendre toutes les informations contenues dans la notice d'utilisation.

Ne pas lire, comprendre et respecter les instructions contenues dans ce document peut endommager le produit, causer des blessures au personnel d'exploitation, ou limiter les performances de l'instrument. Tecan n'est pas responsable des dommages ou des blessures causés par une manipulation incorrecte de l'appareil.

- 3. Respectez tous les avertissements intitulés ATTENTION et PRÉCAUTION de ce document.
- 4. Débranchez toujours l'appareil de l'alimentation principale avant de le nettoyer et de le désinfecter.
- 5. N'ouvrez jamais le boîtier de l'instrument.
- 6. Respectez les mesures de sécurité adaptées du laboratoire, telles que le port de vêtements de protection (par ex. gants, blouse et lunettes de sécurité), et appliquez les procédures de sécurité approuvées du laboratoire.



#### PRÉCAUTION

#### SI LES INSTRUCTIONS DONNÉES DANS CE MANUEL NE SONT PAS SUIVIES CORRECTEMENT, L'INSTRUMENT PEUT ÊTRE ENDOMMAGÉ OU LES PROCÉDURES PEUVENT NE PAS ÊTRE EXÉCUTÉES CORRECTEMENT, ET LA SÉCURITÉ DE L'INSTRUMENT NE PEUT ÊTRE GARANTIE.

On suppose que grâce à leur expérience professionnelle, les opérateurs de l'instrument sont familiarisés avec les précautions de sécurité nécessaires à la manipulation de produits chimiques et biologiques dangereux.

Respectez les législations et règles suivantes :

- Législation nationale sur la sécurité industrielle
- Régulations de prévention des accidents
- Fiches de sécurité des fabricants de réactifs







SELON LES APPLICATIONS, CERTAINES PARTIES DE L'INFINITE F50 PLUS PEUVENT ENTRER EN CONTACT AVEC DES PRODUITS PRÉSENTANT UN RISQUE BIOLOGIQUE OU INFECTIEUX.

ASSUREZ-VOUS QUE SEUL LE PERSONNEL QUALIFIÉ UTILISE L'INSTRUMENT. SI L'INSTRUMENT DOIT ÊTRE RÉPARÉ, DÉPLACÉ OU ÉLIMINÉ, VEILLEZ À TOUJOURS DÉSINFECTER L'INSTRUMENT EN SUIVANT LES INSTRUCTIONS DONNÉES DANS CE DOCUMENT.

RESPECTEZ LES PRÉCAUTIONS DE SÉCURITÉ ADAPTÉES DU LABORATOIRE, TELLES QUE LE PORT DE VÊTEMENTS DE PROTECTION, LORSQUE VOUS TRAVAILLEZ AVEC DES SUBSTANCES POTENTIELLEMENT INFECTIEUSES.

#### AVERTISSEMENT

L'INSTRUMENT RESPECTE LES EXIGENCES EN MATIÈRE D'ÉMISSIONS ET D'IMMUNITÉ DÉTAILLÉES DANS LA NORME IEC 61326-2-6 ; CEPENDANT, L'ENVIRONNEMENT MAGNÉTIQUE DOIT ÊTRE ÉVALUÉ AVANT D'UTILISER L'INSTRUMENT.

IL EST DE LA RESPONSABILITÉ DE L'OPÉRATEUR DE S'ASSURER DU MAINTIEN D'UN ENVIRONNEMENT ÉLECTROMAGNÉTIQUE COMPATIBLE AVEC L'INSTRUMENT, AFIN QUE L'INSTRUMENT FONCTIONNE TEL QUE PRÉVU.

N'UTILISEZ PAS L'INSTRUMENT À PROXIMITÉ DE SOURCES DE RADIATIONS ÉLECTROMAGNÉTIQUES FORTES (PAR EX. SOURCES RF INTENTIONNELLES NON PROTÉGÉES), CAR CELA PEUT GÊNER LE BON FONCTIONNEMENT DE L'INSTRUMENT ET ÉGALEMENT CONDUIRE À DES RÉSULTATS INCORRECTS.





# 2. Général

# 2.1 Utilisation prévue

L'INFINITE F50 PLUS est un lecteur d'absorbance automatisé de microplaques à 96 puits qui inclut le logiciel Magellan. Il est prévu pour une utilisation professionnelle en laboratoire dans les diagnostics in vitro et destiné à mesurer l'absorbance lumineuse (densité optique) de médias liquides homogènes.

L'instrument est prévu pour être utilisé principalement pour réaliser des analyses de diagnostic in vitro sur des échantillons issus du corps humain fournis par des dosages enzymatiques (ELISA). Les informations de diagnostic spécifiques et le type d'échantillon sont définis par le dosage sélectionné.

L'INFINITE F50 PLUS est destiné à la mesure et l'évaluation de dosages qualitatifs, semi-quantitatifs et quantitatifs réalisés conformément aux paramètres de diagnostic programmés et aux spécifications de l'instrument.

Le produit est prévu pour une utilisation professionnelle en laboratoire par du personnel formé. Le produit n'est pas destiné à une utilisation domestique ou par une personne non formée.

Note





Si le lecteur d'absorbance INFINITE F50 PLUS ou le logiciel Magellan sont modifiés de quelque manière, la garantie ne sera plus valide, et l'instrument ne sera plus conforme aux normes en vigueur.

#### Note

L'autorité opérationnelle doit utiliser uniquement des kits de test labellisés CE pour les applications de diagnostic clinique. L'autorité opérationnelle doit s'assurer que la combinaison d'un kit labellisé CE avec le lecteur d'absorbance labellisé INFINITE F50 PLUS et ses options a été conçue pour respecter le règlement IVD 2017/746 ou toute autre régulation nationale ou locale applicable.

Si le lecteur d'absorbance INFINITE F50 PLUS est utilisé pour des applications différentes de celles décrites ci-dessus, ou s'il est utilisé avec un logiciel autre que Magellan, il n'est plus conforme à l'IVD, et l'autorité opérationnelle est responsable de son usage et des validations nécessaires.

# i

Note

Les résultats obtenus en utilisant l'INFINITE F50 PLUS sont déterminés par une utilisation correcte de l'instrument et des microplaques, selon les instructions données dans ce document, ainsi que par les composés liquides utilisés (réactifs, produits chimiques). Les instructions d'utilisation, de stockage et d'application utilisant des échantillons ou des réactifs doivent être strictement respectées. Les résultats doivent donc être interprétés avec prudence.



Note N'ouvrez jamais le boîtier de l'instrument, sous peine de nullité de la garantie.

# 2.2 **Principe de fonctionnement**

L'INFINITE F50 PLUS est un lecteur d'absorbance à 8 canaux destiné à la mesure de la densité optique (DO) des liquides dans des microplaques à 96 puits.

La lumière produite par la source lumineuse DEL traverse un filtre optique dans la roue porte-filtres afin de sélectionner la longueur d'onde nécessaire pour la mesure d'absorbance.

Après avoir traversé le filtre, le rayon lumineux est divisé en huit fibres optiques qui sont focalisées sur huit puits de la microplaque. Les rayons lumineux traversant les puits de la microplaque sont focalisés par huit lentilles optiques puis sont envoyés vers huit photodiodes afin d'être détectés.

Cela signifie que le lecteur INFINITE F50 PLUS mesure huit puits en parallèle, ce qui correspond à une colonne d'une microplaque à 96 puits.

Pour le contrôle et la régulation de l'intensité de lumière de la lampe DEL, une fibre optique supplémentaire (canal de référence) est utilisée, dans laquelle la lumière contourne la microplaque.

Les valeurs de transmission obtenues sont converties en valeurs de densité optique (DO) selon la formule suivante :

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

l<sub>0</sub> = intensité de la lumière incidente

I = intensité de la lumière transmise

La DO est le logarithme de la transmission réciproque.

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



# 2.3 Profils d'utilisateur

# 2.3.1 Utilisateur professionnel – Niveau Administrateur

L'administrateur est une personne disposant d'une formation technique adaptée et des compétences et expériences correspondantes. Si le produit est utilisé comme prévu, la personne est en mesure d'identifier et d'éviter des dangers.

L'administrateur dispose de vastes compétences et est en mesure d'instruire l'utilisateur final ou l'utilisateur routinier dans des protocoles d'essai en rapport avec un produit Tecan dans le cadre de son utilisation prévue.

Des compétences informatiques et de bonnes compétences en anglais sont requises.

### 2.3.2 Utilisateur final ou utilisateur routinier

L'utilisateur final ou l'utilisateur routinier est une personne disposant d'une formation technique adaptée et des compétences et expériences correspondantes. Si le produit est utilisé comme prévu, la personne est en mesure d'identifier et d'éviter des dangers.

Des compétences informatiques et de bonnes compétences linguistiques dans la langue nationale respective sur le site d'installation ou en anglais sont requises.

#### 2.3.3 Techniciens de maintenance

Le technicien de maintenance est une personne disposant d'une formation technique adaptée et des compétences et expériences correspondantes. Si le produit doit être soumis à une révision ou une maintenance, la personne est en mesure d'identifier et d'éviter des dangers.

Des compétences informatiques et de bonnes compétences en anglais sont requises.



Note Les dates de formation, leur durée ainsi que leur fréquence sont disponibles auprès de votre support client.

L'adresse et le numéro de téléphone se trouvent sur Internet : <u>http://www.tecan.com/customersupport</u>



# 3. Prise en main

# 3.1 Déballage et inspection

### 3.1.1 Inspection de l'emballage livré

L'instrument livré inclut :

- Alimentation externe
- Câble d'alimentation
- Câble USB pour connexion à un ordinateur externe
- Notice d'utilisation, fichiers PDF sur support de données
- Clé USB
  - o Logiciel Magellan
  - Notice d'utilisation au format PDF
  - Outils (par ex. Adobe Reader)



Note Pour éviter toute perte de données non désirée ou toute attaque par un virus ou un malware, ne retirez jamais la protection en écriture de la clé USB.



#### PRÉCAUTION

#### LE LECTEUR A ÉTÉ TESTÉ AVEC LE CÂBLE USB FOURNI. SI UN AUTRE CÂBLE USB EST UTILISÉ, IL EST IMPOSSIBLE DE GARANTIR QUE LES PERFORMANCES DE L'INSTRUMENT SERONT CORRECTES.

# 3.1.2 Procédure de déballage

- 1. Vérifiez visuellement que l'emballage n'est pas endommagé avant de l'ouvrir. *Signalez tout dommage immédiatement.*
- Sélectionnez un emplacement où poser l'instrument. L'emplacement doit être plat, sans vibration, à l'abri de la lumière du soleil, et dépourvu de poussière, de solvants et de vapeurs acides. Assurez-vous que la distance entre l'instrument et le mur ou tout autre équipement est d'au moins 5 cm.
- 3. Sortez l'instrument du carton en le soulevant et placez à l'endroit choisi. Soyez précautionneux lorsque vous soulevez l'instrument.
- Procédez à une inspection visuelle de l'instrument, à la recherche de pièces desserrées, tordues ou cassées.
  - Signalez tout dommage immédiatement.
- 5. Comparez le numéro de série de l'instrument, visible sur la plaque située audessous de l'instrument, avec le numéro de série figurant sur l'emballage. *Signalez immédiatement toute incohérence.*
- 6. Vérifiez que les accessoires de l'instrument correspondent à la note accompagnant l'emballage.
- 7. Conservez l'emballage pour pouvoir le réutiliser lors de futurs transports.



# 3.2 Caractéristiques d'alimentation nécessaires

L'instrument détecte automatiquement le voltage fourni. Il n'est donc pas nécessaire de modifier la plage de voltage. Vérifiez les spécifications de voltage et assurez-vous que le voltage fourni à l'instrument correspond aux spécifications suivantes :

Voltage :	
Instrument de base avec adaptateur AC :	100 – 240 V AC, 50/60 HZ
Instrument de base sans adaptateur AC :	24 V DC

Si le voltage suivant n'est pas disponible dans votre pays, veuillez contacter le service client Tecan local.

Ne connectez l'instrument qu'à un système d'alimentation électrique relié à la terre.



PRÉCAUTION N'UTILISEZ PAS L'INSTRUMENT AVEC UNE PLAGE DE VOLTAGE INCORRECTE. SI L'INSTRUMENT EST ALLUMÉ ALORS QU'IL EST CONNECTÉ À UN VOLTAGE INCORRECT, IL SERA ENDOMMAGÉ.



#### PRÉCAUTION

NE REMPLACEZ PAS LES CÂBLES D'ALIMENTATION PRINCIPALE AMOVIBLES PAR DES CÂBLES DE QUALITÉ INSUFFISANTE.

# 3.3 Caractéristiques environnementales nécessaires

L'instrument doit être placé sur une surface plate et égale, dépourvue de poussière, de solvants et de vapeurs acides.

Il est nécessaire d'éviter les vibrations et l'exposition directe à la lumière du soleil pour assurer des résultats corrects.

Température ambiante :	
Fonctionnement	15 °C à 35 °C (59 °F à 95 °F)
Stockage	-30 °C à 60 °C (-22 °F à 140 °F)
Humidité relative :	20% à 80% non condensante à température de fonctionnement



# **3.4 Caractéristiques système nécessaires**

	Minimum	Recommandé
Ordinateur	Ordinateur Windows compatible avec processeur compatible Pentium tournant à 1 GHz (Dual Core)	2 GHz (Dual Core)
Système	Windows 10 (32 bits)	
d'exploitation	Windows 10 (64 bits)	
	Édition : Pro	
Mémoire	Windows 10 (32 bits) : 1 Go RAM	2 Go RAM
	Windows 10 (64 bits) : 2 Go RAM	4 Go RAM
Espace disponible sur le disque dur	3 Go	5 Go
Moniteur	Super VGA Graphics	
Résolution	1024 x 600 et supérieure	1920 x 1080
Profondeur des couleurs	256	
Souris	Souris Microsoft ou dispositif de pointage compatible	
Communication	1 x USB 2.0	2 x USB 2.0 1 x RS232 (Série)
Périnhériques	Windows 10 ·	
i onprioriduoo	Dispositif graphique DirectX 9 avec pilote WDDM 1.0 ou plus récent	
.NET	Microsoft .NET Framework 3.5 :	
	Sous Windows 10, il est demandé à l'utilisateur d'installer la version requise de .NET Framework (version 3.5), si elle n'est pas déjà installée.	
Windows Installer	3.1	
	Si cette version n'est pas installée, le programme d'installation/de mise à niveau l'installera.	
Microsoft Excel	2007, 2010, 2013, 2016 (32 bits)	2010 (32 bits)
	2019 (32 bits)	2019 (32 bits)
	Microsoft Excel 365 (32 bits)	
	Seules les éditions 32 bits sont prises en charge !	
	Les éditions Starter ne sont PAS prises en	
	charge !	

# 3.5 Mettre l'instrument sous tension

Les procédures suivantes décrivent les étapes devant être accomplies avant de mettre l'instrument sous tension.



#### PRÉCAUTION

#### AVANT D'INSTALLER L'INSTRUMENT ET DE LE METTRE SOUS TENSION, IL EST NÉCESSAIRE DE LE LAISSER REPOSER AU MOINS TROIS HEURES DE MANIÈRE À ÉLIMINER TOUT RISQUE DE COURT-CIRCUIT CAUSÉ PAR LA CONDENSATION.

Lorsque les étapes décrites ci-dessus ont été effectuées, l'installation s'effectue selon la procédure suivante :

- 1. Connectez l'instrument à l'ordinateur externe avec le câble USB.
- 2. Assurez-vous que l'interrupteur principal, situé sur le côté gauche de l'instrument, est en position OFF.
- 3. Insérez le câble d'alimentation dans la prise d'alimentation principale du panneau gauche.
- 4. Mettez l'instrument sous tension grâce à l'interrupteur situé sur le panneau gauche.

L'instrument sera prêt à mesurer des microplaques une fois les logiciels installés.



#### Note

Avant de commencer les mesures, assurez-vous que la position A1 de la microplaque est insérée correctement.

Les microplaques ne peuvent être mesurées que sans capot.

Refermez le couvercle du transport de plaque avant de démarrer une mesure, afin d'éviter que la lumière ambiante n'influe sur les résultats.



Note Toujours décharger la microplaque du lecteur immédiatement après la fin de la mesure.



# 3.6 Logiciel

# 3.6.1 Introduction/ Utilisation prévue

Le logiciel de contrôle d'instrument et d'analyse des données **Magellan** est livré avec l'instrument.

Magellan est un **logiciel de contrôle de lecteur et d'analyse de données** universel, destiné à analyser les données générées par des tests de microplaques utilisant des appareils de mesure Tecan.

Magellan est disponible en deux versions :

- Magellan et
- Magellan Tracker

Le logiciel **Magellan** est prévu pour effectuer des mesures de point final, cinétiques et multi-étiquetages avec l'instrument INFINITE F50 PLUS, conformément à son utilisation prévue ; voir le chapitre 2.1 Utilisation prévue.

**Magellan Tracker** offre toutes les fonctionnalités nécessaires pour entrer en conformité avec la régulation 21 CFR partie 11 de la FDA, en plus de la fonctionnalité de Magellan.



Note

Veuillez noter qu'une installation correcte de l'instrument et du logiciel Magellan seule ne garantit pas d'être en conformité avec les régulations. Des règles adaptées aux processus et aux procédures opérationnelles permanentes, notamment de validation et de contrôle qualité, doivent également être mises en place.

# 3.6.2 Procédure d'installation

Pour installer le logiciel, insérez la clé USB dans le port USB, et procédez comme suit :

- L'Assistant d'installation de Magellan devrait démarrer automatiquement et vous guider tout au long du processus d'installation. Si ce n'est pas le cas, veuillez lancer le fichier "E:\Tecan.exe" (si E est la lettre attribuée à la clé USB).
- Sélectionnez 'Magellan software' et en fonction de la version que vous avez commandée – 'Install Magellan' ou 'Install Magellan Tracker' pour lancer la procédure d'installation, puis suivez l'assistant.
- 3. Cliquez sur Installer pour lancer la procédure d'installation du logiciel.
- 4. Cliquez sur **I accept** the terms of the license agreement puis sur **Next** pour continuer.
- 5. La page **Informations client** apparaît alors : veuillez entrer votre nom et celui de votre organisation.
- 6. La Page de configuration apparaît : choisissez votre langue.
- 7. Page : Use for regulated environments : cliquez sur Next pour continuer.
- 8. Cliquez sur Installer pour lancer l'installation.
- 9. Cliquez sur **Terminer** pour finir l'installation et fermer le programme d'installation.

Le logiciel peut être lancé via le menu **Démarrer** de Windows, en sélectionnant **Magellan** dans le groupe de programmes **Tecan**.



Note Il est très important que la personne qui installe le logiciel dispose de privilèges d'administrateurs sur l'ordinateur.





Note Magellan V7.5 ne peut pas être installé parallèlement à d'autres versions du logiciel Magellan.

Par défaut, tous les types de fichiers associés à Magellan sont stockés dans des sous-dossiers correspondants du répertoire suivant :

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

## 3.6.3 Qualification de l'installation

Vérifiez que l'installation de Magellan a réussi grâce au programme automatique de qualification de l'installation :

Lancez *TecanIQ.exe* depuis le chemin d'installation par défaut (C: *\Program Files\TecanWagellan*) ou depuis le **menu Démarrer Windows : Démarrer > Programmes > Tecan > MagellanIQ.** 

Cliquez sur **Vérifier** pour lancer la qualification de l'installation. Tous les composants installés devraient avoir le statut **OK**. Veuillez contacter votre représentant Tecan local si un problème potentiel est signalé.

Pour fermer le programme de qualification de l'installation, cliquez sur **Annuler** ou sur **Quitter** 



Note

La qualification de l'installation doit être répétée à chaque installation ou mise à jour vers une nouvelle version de Magellan.

# 3.6.4 Commencer à travailler avec Magellan

Le type d'interface principal de Magellan est l'assistant.

Les assistants de Magellan représentent les modules de flux de travail, qui sont des guides étape par étape pour effectuer des procédures complexes.

Parfois, des menus sont disponibles dans la barre d'en-tête. Le **Menu** offre un moyen conventionnel d'utiliser le logiciel : l'élément de menu approprié est sélectionné dans les menus principaux. Toutes les actions en découlant sont démarrées instantanément, ou une boîte de dialogue est affichée, dans laquelle il est possible d'effectuer des sélections supplémentaires ou des saisies.



#### Interface utilisateur – liste des assistants

Après le démarrage de Magellan, la Liste des assistants apparaît :

Liste des assistants - Magellan Tracker		
0	Que voulez-vous faire?	
Démarrer mesure Évaluer les résultats	<ul> <li>L'assistant Démarrer la mesure vous aide à effectuer une mesure. Vous pouvez utiliser une méthode ou obtenir des données brutes</li> </ul>	
Attacher une signature Créer/éditer une liste ID	•	
d'echantilion Créer/éditer une méthode	•	
Quitter Magellan		

Chaque assistant peut être démarré par un double clic ou en le sélectionnant et

en cliquant sur le bouton Suivant

#### Assistant Démarrer la mesure

L'assistant Démarrer la mesure comprend les options suivantes :

- Obtenir les données brutes est utilisé pour générer rapidement et facilement des données brutes en définissant les paramètres de mesure requis et en démarrant une mesure.
- Utiliser méthode prédéfinie est utilisé pour effectuer des mesures basées sur des méthodes précédemment définies.
- Démarrer favoris est utilisé pour sélectionner l'une des méthodes les plus fréquemment utilisées dans la liste d'icônes numérotées.

Une fois la mesure terminée, un fichier espace de travail est créé.

#### Assistant Évaluer les résultats

L'assistant **Évaluer les résultats** est utilisé pour visualiser les données brutes et pour évaluer les résultats. Les paramètres d'évaluation peuvent être visualisés et les données réévaluées.

#### Assistant Attacher une signature

L'assistant **Attacher une signature** est utilisé pour signer les fichiers de méthode et d'espace de travail. Cette fonctionnalité n'est disponible qu'avec Magellan Tracker.

#### Assistant Créer/éditer une liste ID d'échantillon

L'assistant **Créer/éditer une liste ID d'échantillon** est utilisé pour créer une nouvelle liste ID d'échantillon et éditer des listes existantes.

#### Assistant Créer/éditer une méthode

L'assistant **Créer/éditer une méthode** est utilisé pour définir ou éditer des méthodes.

#### 3. Prise en main





#### Note Pour des informations détaillées sur le logiciel, veuillez consulter la notice d'utilisation de Magellan.

Veuillez noter que certaines fonctionnalités décrites dans la notice d'utilisation de Magellan peuvent ne pas concerner (être indisponibles dans) Magellan V7.5 en combinaison avec l'INFINITE F50 Plus. Cependant, toutes les informations nécessaires sont présentes dans cette notice d'utilisation.



Note Vous trouverez un exemple détaillé de mesure ELISA au chapitre 6 Exemple d'application.

Les fichiers de méthode Sunrise créés avec Magellan V7.x ou supérieur peuvent être ouverts avec Magellan V7.5; les paramètres de mesure sont convertis automatiquement.

Les méthodes Sunrise créées avec des versions plus anciennes doivent être converties en utilisant l'option "Convertir à partir de" du menu Divers/Gestion des fichiers.



# 3.7 Magellan - Éditeur de paramètres de mesure

L'assistant **Editeur de paramètres de mesure** est utilisé pour mettre en place des flux de travail. Chaque flux de travail est créé facilement en glissant et déposant les étapes du processus sous forme de séquence, selon l'application. Le flux de travail de l'application peut ensuite être visualisé par l'utilisateur dans le volet flux de travail. Chaque étape du processus (élément de programme) peut être copiée et collée (en utilisant les raccourcis standards de Windows, **Ctrl-C**, **Ctrl-V** ou le menu contextuel de la souris), et déplacée à l'endroit souhaité dans le flux de travail.



L'Éditeur de paramètres de mesure comprend les éléments suivants, qui sont décrits en détail dans les prochains chapitres :

- Barre de contrôle
- Volet Flux de travail
- Volet d'info



### 3.7.1 Barre de contrôle

La **Barre de contrôle** est divisée en cinq sections. Chaque section contient des éléments de programme utilisés pour créer un flux de travail individuel.

Pour créer un flux de travail, double cliquez sur l'élément de programme sélectionné, ou glissez-le et déposez-le dans le volet Flux de travail.

Les éléments de programme suivants sont disponibles avec l'INFINITE F50 PLUS :

Outils Iaboratoire	Plaque Partie de la plaque
Mesures	Absorbance
Actions	Agitation Déplacer plaque
Cinétique	Cycle cinétique Condition cinétique
Divers	Commentaire Requête de l'utilisateur Attendre (temps) Incubation

#### **Outils laboratoire**

#### Plaque

L'élément de programme **Plaque** est utilisé pour sélectionner un format de plaque dans la liste déroulante **Définition de plaque**. Cliquez sur **Détails**... pour visualiser plus d'informations sur la plaque sélectionnée.

🔷 🔻 Plaque		1
Définition de la plaque :	[COS96ft] - Corning 96 Flat Transparent	V Détails
		<u>Utiliser une partie de la plaque</u>

#### Partie de la plaque

Par défaut, l'élément de programme **Partie de la plaque** est réduit. Lorsqu'il est agrandi en cliquant sur **I**, une microplaque à 96 puits s'affiche. Pour mesurer des puits individuels, cliquez sur le puits choisi ; pour mesurer un intervalle de puits, glissez un cadre autour de l'intervalle choisi. Cliquer sur **Détails** permet de zoomer sur l'aperçu de la plaque.

췤 🔻 Partie de la plaque		2
1       2       3       4       5       6       7       8       9       10       11       12         A       • <th>Détails</th> <th></th>	Détails	



#### Parties de plaque séparées

Des parties plaque séparées peuvent être sélectionnées :



Il est possible de sélectionner un second intervalle de puits en appuyant sur la touche **Contrôle** du clavier et en glissant un cadre sur les puits à sélectionner.

#### Mesures

#### Absorbance

L'élément de programme **Absorbance** est utilisé pour effectuer des mesures d'absorbance. Entrez ou sélectionnez les paramètres respectifs.

Deux listes déroulantes contiennent les longueurs d'onde des filtres de mesure et de référence disponibles, en fonction des filtres d'absorbance insérés. Si les listes déroulantes sont vides, les filtres n'ont pas été définis.

🖕 🔻 Absorba	ance				3
-Longueur d'onde	)	Étiquet	tte		
Mesure:	492 nm 🗸 🗸	Nom:	Étiquette1	*	
Référence:	405 nm 💌				

#### Actions

#### Agitation

Sélectionnez l'élément de programme **Agitation** si la plaque doit être agitée, soit avant la mesure, soit entre les cycles cinétiques.

🧙 🔻 Agitation		4
Paramètre Durée: 1 📚 sec Inte	nsité: Wide Amplitude: 14,1 mm Fréquence: 2,1 Hz	Attendre guelgues secondes

Entrez les paramètres respectifs :

Durée	Entrez la durée du processus d'agitation.
Intensité	Entrez le mode d'agitation choisi. L'amplitude et la fréquence sont affichées lors du choix d'un mode d'agitation.

Consultez le chapitre 4.1.1 Agitation de la microplaque pour les modes d'agitation disponibles.



#### 3. Prise en main

Cliquer sur le lien <u>Attendre quelques secondes</u> insère un nouvel élément de programme. Voir la page 25 pour plus de détails.

#### Cinétique

#### Cycle cinétique

Utilisez l'élément de programme **Cycle cinétique** pour effectuer plusieurs mesures consécutives, qui peuvent être exécutées à des intervalles donnés.

🎧 ▼ Cycle cinétique		3
Cycles ···· 2 2 • Nombre de cycles : 2 2 Durée	Intervale cinétique         ✓ Utiliser intervale cinétique:         ④ Temps :       00:01:00 ♀) (hh:mm:ss)         ● Temps :       60000 ♀) ms	

Entrez les paramètres respectifs :

Cycles	<ul> <li>Nombre de cycles : Entrez un nombre ou cliquez sur les flèches haut ou bas pour définir le nombre d'étapes de mesure (2 – 1000 cycles).</li> <li>Durée : Entrée la durée au format hh:mm:ss</li> </ul>
Intervalle cinétique	<b>Utiliser intervalle cinétique</b> : Entrez l'intervalle temporel (hh:mm:ss ou ms).

#### **Condition cinétique**

Utilisez l'élément de programme **Condition cinétique** pour définir quelles actions doivent être exécutées lors d'un certain cycle.

🙊 🔻 Condition cinétique	4
Condition Exécuter les commandes au cycle : 2	

Si 2 est entré pour **Exécuter commande au cycle** pour une mesure cinétique comprenant, par exemple, une étape **Agitation**, l'agitation sera effectuée au cycle 2.



Note

Les conditions cinétiques telles que l'agitation doivent être insérées immédiatement après un élément de programme Cycle cinétique, afin d'assurer une répétabilité maximale des résultats. Il est conseillé aux utilisateurs de mettre en place des scripts adaptés avant de faire les mesures, et d'utiliser le même script pour toutes les mesures cinétiques similaires afin d'obtenir des résultats comparables.

**Divers** 

#### Commentaire

Utilisez l'élément de programme **Commentaire** pour entrer une remarque ou un commentaire concernant la mesure en cours dans le champ de texte.

💷 🔻 Commentaire	6
Commentaire :	



#### Requête de l'utilisateur

L'élément de programme **Requête de l'utilisateur** informe l'opérateur de l'instrument qu'il doit exécuter une action définie à un certain moment du flux de travail.

🚯 🔻 Requête de l'utilisateur	7
Texte :	

Si, par exemple, l'élément de programme **Déplacer plaque** est utilisé pour sortir la plaque afin d'effectuer une action, le texte entré doit informer l'opérateur qu'il doit effectuer cette action. Une boîte de dialogue affiche le message, et le processus de mesure s'arrête jusqu'à ce qu'**OK** soit cliqué.

#### Attendre (horloge)

Utilisez l'élément de programme **Attendre (horloge)** pour définir une période d'attente avant l'exécution de l'étape suivante du flux de travail.

#### Entrez la durée choisie dans le champ Attente.

🮯 🔻 Attendre (horloge)		5
Attente : 00:01:00 🗘 (hh:mm:ss)	Options Attendre l'injection Ignorer l'attente pour le dernier cycle cinétique	

Entrez les paramètres respectifs :

Horloge	Entrez l'Attente (hh:mm:ss)
Options	Ignorer l'attente pour le dernier cycle cinétique : Si l'étape du programme Attendre (horloge) est la dernière action d'une exécution cinétique, le temps d'attente sera ignoré lors du dernier cycle.

#### Incubation

🥝 ▼ Incubation	4
Horloge Temps d'incubation: 00:01:00 📚 (hh:mm:ss)	

Entrez les paramètres appropriés pour l'incubation :

Temps	Entrez le temps total (min. 5s)
d'incubation	



# 3.7.2 Volet Flux de travail

Le **Volet flux de travail** est la fenêtre où le script de mesure est affiché et où les paramètres sont définis et édités.

Il y a deux manières d'insérer un élément de programme de la **Barre de contrôle** dans le **Volet Flux de travail** :

- Sélectionnez un élément de programme dans la Barre de contrôle ; si vous double cliquez dessus, il est inséré dans le Volet Flux de travail directement après l'élément de programme précédent.
- Cliquez sur un élément de programme dans la **Barre de contrôle** et glissez-le dans le **Volet Flux de travail** jusqu'à la position souhaitée.

Les éléments de programme sont numérotés en fonction de leur position dans la séquence.

Une fois qu'un élément de programme a été inséré dans le **Volet Flux de travail**, les réglages et les paramètres de cet élément peuvent être entrés ou édités.

Les éléments de programme contenus dans le **Volet Flux de travail** peuvent être réduits pour ne montrer que les informations les plus importantes, ou agrandis pour permettre d'accéder à toutes les fonctions éditables. Cliquez sur l'un des triangles situés à côté de l'élément de programme, vou , pour passer d'un mode d'affichage à l'autre.

Par défaut, l'éditeur de paramètres de mesure se lance avec l'élément **Plaque**, l'élément **Partie de plaque** (réduit), et un élément **Absorbance** dans le **Volet Flux de travail**.

Les éléments de programme actuellement sélectionnés dans le **Volet Flux de travail** s'affichent avec une ligne jaune le long de leur bordure supérieure.

Si un élément de programme contient des erreurs ou n'est pas valide dans le flux de travail actuel, il sera signalé par une marque d'erreur, et son numéro d'élément sera surligné en rouge. Des informations sur l'erreur s'affichent dans le **Volet d'info**. Si le flux de travail contient des erreurs, les paramètres de mesure ne peuvent pas être choisis.

#### Hiérarchie des éléments

La hiérarchie des éléments dans le Volet Flux de travail est la suivante :

- 1. Plaque
- 2. Partie de la plaque (Intervalle)

Toute étape de mesure souhaitée peut être insérée directement après un élément de plaque ou d'intervalle. Utilisez **Libérer** et **Indenter** pour modifier la séquence d'exécution d'un composant à une seule bande. Sélectionnez un élément dans le **Volet Flux de travail**, cliquez avec le bouton droit de la souris, et sélectionnez **Libérer** ou **Indenter**.

D'autres éléments de la **Barre de contrôle** peuvent être insérés dans la hiérarchie d'un flux de travail :

Le premier élément **Intervalle** est inséré directement après l'élément **Plaque** ; tous les éléments **Intervalle** suivants peuvent ensuite être insérés.

Les étapes cinétiques sont possibles avec un élément Plaque ou Intervalle.

Des étapes **Requête utilisateur**, **Commentaire** et **Attendre** sont possibles dans un élément **Plaque** ou **Intervalle**.

#### 3.7.3 Volet d'info

Le **Volet d'info** situé à droite de l'écran affiche les informations concernant l'élément de programme actuellement sélectionné. Tout avertissement ou erreur y est affiché.



# 3.8 Magellan - Définir les mesures

Le chapitre qui suit présente des exemples illustrant la définition de différentes mesures.

#### 3.8.1 Définir les mesures de point final

L'exemple qui suit décrit une **Mesure de point final d'absorbance** dans tous les puits d'une microplaque à 96 puits.

- 1. Sélectionnez une microplaque à 96 puits dans la liste déroulante **Définition de plaque**.
- 2. Par défaut, tous les puits d'une microplaque à 96 puits sont sélectionnés pour la mesure.
- 3. Entrez la mesure choisie et les longueurs d'onde de référence.

🔷 ▼ Plaque	1
Définition de la plaque : [COS96ft] - Corning 96 Flat Transparent 🕑 Détails	
Utiliser une partie de la pla	que
▼ Partie de la plaque         1       2       3       4       5       6       7       8       9       10       11       10         A       ● <t< td=""><td>2</td></t<>	2
H Absorbance	3
Référence: 620 nm	



# 3.8.2 Définir des mesures multi-étiquetage

L'exemple qui suit décrit une **Mesure d'absorbance multi-étiquetage** dans un intervalle défini d'une microplaque à 96 puits (A1:E7). Trois étiquettes d'absorbance seront mesurées.

- 1. Sélectionnez une microplaque à 96 puits dans la liste déroulante **Définition de plaque**.
- 2. Par défaut, tous les puits d'une microplaque à 96 puits sont sélectionnés pour la mesure.

Cliquez sur pour agrandir l'élément **Partie de plaque**. Ensuite, sélectionnez l'intervalle désiré (A1:E7).

- 3. Entrez la longueur d'onde de mesure souhaitée.
- 4. Insérez 2 éléments **Absorbance** de plus et entrez les longueurs d'onde de mesure.

🔶 🔻 Plaque			1
Définition de la plaque	e : [COS96ft] - Corning 96 Flat Transparent	▼	Détails
	L	Utiliser une parti	e de la plaque
🎪 🔻 Pa	rtie de la plaque		2
1 2 B 0 C 0 E 0 G 0 H	3 4 5 6 7 8 9 10 11 12 • • • • • • • • • • • • • • • • • • •	is	
	▼ Absorbance		3
Long	gueur d'onde	Étiquette	
Mes	ure: 405 nm 💌	Nom: Étiquette1	
R	Reference: 405 nm		
	▼ Absorbance		4
Long	gueur d'onde	-Étiquette	
Mes	ure: 492 nm 💌	Nom: Étiquette2	
R	Référence: 405 nm		
	▼ Absorbance		5
Long	gueur d'onde	Étiquette	
Mes	sure: 620 nm 💌	Nom: Étiquette3	
R	Référence: 405 nm		



# 3.8.3 Définir des mesures cinétiques

L'exemple qui suit décrit une mesure cinétique d'une microplaque à 96 puits.

- 1. Sélectionnez une microplaque à 96 puits dans la liste déroulante **Définition de plaque**.
- 2. Insérez un élément de programme **Cycle cinétique** entre la partie de plaque et l'élément absorbance.
- 3. Cycles/Nombre de cycles : 50
- 4. Intervalle cinétique (intervalle entre les mesures) : sélectionnez **Utiliser** intervalle cinétique et entrez : 2 minutes 30 secondes.
- 5. Définissez l'élément **Absorbance** en entrant la longueur d'onde de mesure souhaitée.

🔶 🔻 Plaque		1
Définition de la plaque :	[COS96ft] - Corning 96 Flat Transparent	Détails Utiliser une partie de la plaque
1         2         3           A         -         3           B         -         6           C         -         6           D         -         6           F         -         6           H         -         6	de la plaque	2
Gov ← C Cycles ⊙ Nombre ○ Durée	Eycle cinétique a de cycles : 50 ♥ Utiliser intervalle cinétique : ④ Temps : 00:02:30 ♥ (hh:mm:ss) ◯ Temps : 150000 ♥ ms	3
	▼ Absorbance ongueur d'onde esure: 192 nm ▼ Référence: 105 nm ▼	4
		05



### 3.8.4 Indentation et libération des éléments de programmes

Choisir d'indenter/libérer un élément de programme modifiera le flux de travail de l'instrument pendant les mesures.

Les actions de tous les éléments de programme ayant la même indentation sont effectuées consécutivement. La seule dépendance entre ces éléments de programme est que l'action suivante démarre directement après que la précédente s'est terminée.

Un élément de programme plus indenté que le précédent élément de programme signale une dépendance entre les deux éléments de programme. Cela signifie que les paramètres définis dans le premier élément de programme sont également actifs pour le second élément de programme (indenté).

Voici un exemple montrant comment définir une **Cinétique multi-étiquetage** avec deux **Etiquettes d'absorbance**. L'exemple montre que les deux **éléments de programme** Absorbance dépendent de l'élément de programme **Cycle cinétique**, qui dépend de l'**élément de programme** partie de la plaque, qui dépend de l'élément de programme **Plaque**. Définissez les paramètres comme suit pour cet exemple :

- 1. Plaque : par ex. Greiner 96 transparente à fond plat
- 2. Cycle cinétique/Nombre de cycles : 5
- 3. Absorbance/longueur d'onde étiquette 1 : 450 nm
- 4. Absorbance/longueur d'onde étiquette 2 : 492 nm

Le Volet Flux de travail s'affiche comme sur cette capture d'écran :

🔷 🔻 Plaque			1
Définition de la plaque : [GRE96ft]	Greiner 96 Flat Transparent		V Détails
			Utiliser une partie de la plaque
🔹 🔻 Partie de la plag	ue		2
1 2 3 4 5 6 A C C C C C C C C C C	7 8 9 10 11 12 Détails.		
🎧 🔻 Cycle cinét	ique		3
Ovdes   Nombre de cycles :  Durée	5 ¢	ntervalle cinétique ] Utiliser intervalle cinétique	
👌 🔻 Abso	rbance		4
-Longueur d'or	nde	Étiquette	
Mesure:	450 nm 💌	Nom: Étiquette1	
Abso	rbance		5
-Longueur d'or	nde	Étiquette	
Mesure:	492 nm 💌	Nom: Étiquette2	
Référence	405 nm 💌		
			203

La définition ci-dessus aboutit au flux de travail suivant :

L'absorbance de tous les puits de la microplaque à 96 puits est d'abord mesurée à 450 nm, puis à 492 nm. Les deux mesures d'absorbance sont effectuées pendant 5 cycles cinétiques.



Libérer le second élément de programme **Absorbance**, de manière à l'aligner avec l'élément **Cycle cinétique**, modifie le flux de travail. Sélectionnez le second élément de programme **Absorbance** et cliquez du bouton droit de la souris. Sélectionnez **Libérer la bande** dans le menu contextuel. Le **Volet Flux de travail** s'affiche comme sur cette capture d'écran :

🔷 🔻 Plaque	1
Définition de la plaque : [GRE96ft] - Greiner 96 Flat Transparent	Détails le la plaque
i → Partie de la plaque	2
1       2       3       4       5       6       7       8       9       10       11       12         B       Image: Constraint of the state	
Q ▼ Cycle cinétique	3
Cycles       Intervale cinétique         Nombre de cycles :       5 \$         Durée       Utiliser intervale cinétique	
borbance	4
Longueur d'onde       Mesure:       450 nm       Référence:       405 nm	
Absorbance	5
Longueur d'onde Mesure: 492 nm  Référence: 405 nm	

Dans ce flux de travail, une mesure **Cinétique d'absorbance** avec 5 cycles est effectuée pour la première absorbance à 450 nm ; à la fin de cette boucle, la mesure de **Point final d'absorbance** à 492 nm a lieu.



# 3.9 **Optimisation des performances**

L'instrument a été entièrement testé en usine pour assurer que ses performances se situent dans les limites spécifiées (voir 4.4.2 Spécifications des mesures pour plus de détails).

Respecter les recommandations ci-dessous assure une précision maximale à l'instrument.

#### 3.9.1 Emplacement de l'instrument

L'instrument doit être positionné de manière appropriée (voir le chapitre 3.3 Caractéristiques environnementales nécessaires pour plus de détails).

# 3.9.2 Procédure opérationnelle

#### Généralités

- Il est recommandé de suivre les procédures opérationnelles permanentes des dosages effectués.
- Pour obtenir la meilleure répétabilité possible, la longueur d'onde de mesure doit correspondre à la longueur d'onde d'absorbance maximum de la solution étudiée.

Il est important d'utiliser la longueur d'onde d'absorbance maximum si la courbe d'absorbance de l'échantillon est au-dessus d'une bande de longueur d'onde étroite.

Sachez que les mesures effectuées sur la pente d'un pic d'absorbance limiteront la précision des valeurs de DO.

- Après avoir mesuré chaque microplaque, référez-vous au kit de test pour obtenir des informations sur la procédure de validation.
- Utilisez les filtres d'absorbance recommandés pour l'INFINITE F50 PLUS.

#### Microplaques

 L'instrument peut être utilisé avec les types de microplaques présentés au chapitre 4.4.3 Microplaques. Les résultats sont meilleurs en utilisant des microplaques à fond plat. Les résultats des mesures peuvent varier en fonction du type de microplaque utilisé.

Soyez particulièrement prudent lors de l'utilisation de microplaques à fond en C, U ou V ou de plaques à barrettes, car il est possible que les résultats des mesures diffèrent des spécifications décrites dans ce document.

Assurez-vous que le type de microplaque utilisé avec le lecteur d'absorbance INFINITE F50 PLUS convient à l'application qui en est faite.

- N'utilisez que des microplaques parfaitement propres.
- Ne laissez pas de poussière se déposer sur les solutions ou sur la microplaque pendant la période d'incubation préalable à la mesure.
   Il est recommandé d'utiliser un couvercle de protection chaque fois qu'une microplaque est sortie de l'instrument.
- Les inexactitudes dans la quantité de solution pipettée ont un effet important sur les résultats obtenus, lorsque de faibles quantités de solution sont utilisées.
- La forme du ménisque de la solution peut créer des inexactitudes dans les résultats, en particulier lorsque de faibles quantités de solution sont utilisées.



# 4. Fonctionnalités de l'instrument

# 4.1 Fonctionnalités de l'instrument

Les modes de mesure d'absorbance suivants sont disponibles sur l'INFINITE F50 PLUS : mesures de point final, cinétiques, et multi-étiquetages.

### 4.1.1 Agitation de la microplaque

L'INFINITE F50 PLUS est capable d'agiter la microplaque avant la mesure. La microplaque peut également être agitée entre chaque cycle de mesure cinétique. Utilisez Magellan pour définir le mode d'agitation.



#### PRÉCAUTION LORS DE L'AGITATION DE MICROPLAQUE, IL EST POSSIBLE QUE LES PUITS VERSENT S'ILS SONT REMPLIS AVEC UN VOLUME TROP IMPORTANT.

Modes d'agitation de l'INFINITE F50 PLUS :

Mode d'agitation	Amplitude d'agitation	Fréquence d'agitation
ÉLEVÉ	2,8 mm	12,3 Hz
NORMAL	4,4 mm	9,2 Hz
FAIBLE	4,4 mm	7,8 Hz
AMPLITUDE	14,1 mm	2,0 Hz

# 4.2 Description de l'instrument

L'illustration ci-dessous montre les composants de l'instrument.



Position du puits A1 Transport de plaque



#### 4. Fonctionnalités de l'instrument

La DEL d'état donne des informations sur le statut de l'instrument :

- Vert clignotant : l'instrument n'est pas connecté à Magellan
- Vert : l'instrument est connecté et prêt à mesurer
- Rouge : une mesure est en cours

Sur le panneau gauche de l'instrument se trouvent le port USB, l'interrupteur principal et la prise d'alimentation principale.



La plaque d'identification est fixée sous l'instrument.

Exemple de plaque d'identification



Le contenu de la plaque d'identification (nom de modèle et code de l'article) peut changer en fonction du modèle spécifique.

Pour une vue d'ensemble des divers instruments pour lesquels ces instructions d'utilisation s'appliquent, voir la Déclaration de Conformité (Declaration of Conformity) à la dernière page de document.



# 4.3 Description de la roue porte-filtres

La roue porte-filtres standard de l'INFINITE F50 PLUS est livrée avec quatre filtres d'interférences à bande étroite de longueur d'onde fixe (405, 450, 620 et 492 nm). Il est possible d'équiper la roue d'un total de 8 filtres. Veuillez contacter votre représentant Tecan local pour connaître les filtres supplémentaires disponibles.

Les filtres de la roue porte-filtres standard sont montés comme présenté cidessous :

Position du filtre	Longueur d'onde du filtre
1	405 nm
2	450 nm
3	620 nm
4	492 nm
5 - 8	Emplacements vides pour filtres



Lorsqu'une longueur d'onde est sélectionnée pour une mesure, le filtre correspondant est amené vers le rayon lumineux grâce à un déplacement de la roue porte-filtres vers la position correspondante.



Note Pour plus d'informations sur la définition d'un nouveau filtre, voir 7.5.2 Définir les filtres.



# 4.4 Spécifications de l'instrument

Les tableaux ci-dessous contiennent les spécifications du lecteur d'absorbance INFINITE F50 PLUS.

# 4.4.1 Spécifications générales

PARAMETRES	CARACTERISTIQUES
Entrée d'alimentation principale Alimentation externe	Alimentation : Instrument de base avec adaptateur AC : 100 - 240 V AC, 50/60 Hz, max. 1,2 A (détection automatique, survoltage catégorie II) Instrument de base sans adaptateur AC : 24 V DC (survoltage catégorie I)
Consommation INFINITE F50 PLUS	Mode veille : env. 12 VA Mode opérationnel : max. 30 VA
Dimensions extérieures	Largeur :         34,7 cm (13,66 pouces)           Profondeur :         18,9 cm (7,44 pouces)           Hauteur :         13,4 cm (5,28 pouces)
Poids	2,6 kg (alimentation incluse)
Température ambiante :	
Fonctionnement	15 °C à 35 °C (59 °F à 95 °F)
Stockage	-30 °C à 60 °C (-22 °F à 140 °F)
Humidité relative	20 % à 80 %
Degré de pollution	2
Méthode d'élimination des déchets	Déchets contaminés
Environnement	Voir le chapitre 3.3 Caractéristiques environnementales nécessaires pour plus d'informations.


# 4.4.2 Spécifications des mesures

PARAMETRES	CARACTERISTIQUES
Temps de mesure : longueur d'onde simple longueur d'onde double	< 15 secondes < 20 secondes
Intervalle de longueur d'onde : Standard	400 - 750 nm
Intervalle de mesure : 400 - 750 nm	0 - 4,000 DO
Résolution :	0,0001 DO
Exactitude : 450, 492 nm 0,000 - 2,000 DO 2,000 - 3,000 DO	≤ (1.0 % + 0.010 OD)* ≤ (1.5 % + 0.010 OD)*
Précision : 450, 492 nm 0,000 - 2,000 DO 2,000 - 3,000 DO	≤ (0.5 % + 0.005 OD)* ≤ (1.0 % + 0.005 OD)*
Linéarité : 450, 492 nm 0,000 - 2,000 DO 2,000 - 3,000 DO	≤ 1 % ≤ 1.5 %
Sélection de longueur d'onde : Filtre standard	Filtres d'interférence de bande- passante étroite. Une roue porte- filtres peut accueillir jusqu'à 8 filtres.
Précision de la longueur d'onde du filtre :	Longueur d'onde centrale ± 2 nm
Longueur d'onde du filtre : A 50% de transmission	10 ± 2 nm
Source lumineuse :	DEL
Interface informatique :	USB

Tous les appareils connectés doivent être approuvés et listés par la norme IEC 60950-1 Matériels de traitement de l'information – Sécurité et les standards locaux équivalents.

\* supérieure ou égale à x % de la valeur de mesure plus la valeur de DO correspondante



## 4.4.3 Microplaques

Toutes les microplaques à 96 puits et à fond transparent (plat, en C, U, et V ; y compris les microplaques à barrettes) conformes aux standards suivants peuvent être utilisées avec le lecteur d'absorbance INFINITE F50 PLUS :

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



#### PRÉCAUTION N'UTILISEZ QUE DES MICROPLAQUES SANS COUVERCLE, ET N'UTILISEZ PAS DE MICROPLAQUES D'UNE HAUTEUR SUPÉRIEURE À 15,2 MM.

#### Manipulation de la microplaque

N'insérez ou ne retirez la microplaque que lorsque le transport de plaque est entièrement éjecté (voir l'illustration ci-dessous) et que le moteur du transport de plaque est inactif. N'ouvrez pas le capot du boîtier lorsque la DEL d'état est allumée en rouge.



#### AVERTISSEMENT PORTEZ TOUJOURS DES GANTS À USAGE UNIQUE ET DES VÊTEMENTS DE PROTECTION LORSQUE VOUS MANIPULEZ LA MICROPLAQUE.



Position du puits A1 Transport de plaque – entièrement éjecté

# 4.5 Accessoires de l'instrument

La liste ci-dessous présente les accessoires disponibles en option pour l'INFINITE F50 PLUS, qui peuvent être commandés en supplément :

- Filtres supplémentaires
- Outil de montage de filtre
- Plaque MultiCheck<sup>™</sup> pour gamme INFINITE F50

Pour plus d'informations et connaître la disponibilité dans votre pays, veuillez contacter votre représentant Tecan local.



# 5. Contrôle qualité

# 5.1 Introduction



#### PRÉCAUTION SI LES PERFORMANCES ANALYTIQUES DE L'INFINITE F50 PLUS SONT MISES EN DOUTE, SUIVEZ LES INSTRUCTIONS FOURNIES POUR LE CONTRÔLE QUALITÉ OU CONTACTEZ LE CENTRE D'ASSISTANCE TECAN LE PLUS PROCHE.

Ce chapitre fournit des informations sur la procédure d'auto-vérification de l'instrument et des instructions permettant de vérifier facilement la qualité opérationnelle.

# 5.2 Procédure d'auto-vérification

Lorsque l'INFINITE F50 PLUS est connecté au logiciel de contrôle de lecteur Magellan, les moteurs et les senseurs sont vérifiés, et le transport de plaque et la roue porte-filtres sont initialisés.

Avant chaque mesure, une procédure d'auto-vérification de calibrage est effectuée, afin de s'assurer que l'instrument fonctionne correctement et de calibrer le système optique.

# 5.3 Qualification opérationnelle (QO)

Les tests suivants peuvent être effectués afin de s'assurer que l'instrument fonctionne correctement et que les résultats obtenus sont exacts.

La répétabilité et l'exactitude de l'instrument peuvent varier selon le type de solution et de microplaque utilisé.

Pour éliminer cet effet, les instruments sont testés en usine avec une plaque de calibrage, qui supprime l'influence de la solution et toute variation due à la position de la microplaque lors de la mesure.

### 5.3.1 Test MultiCheck

Le test MultiCheck permet de vérifier automatiquement les performances du lecteur, notamment l'exactitude, la linéarité, la précision, et l'alignement avec les standards traçables NIST.

### 5.3.2 Test de microplaque

Si les densités optiques des puits de la microplaque ne sont pas cohérentes, les résultats obtenus avec ce type de microplaque seront influencés.

Cette incohérence peut être vérifiée en lisant une microplaque vide.

Les valeurs de DO obtenues par la mesure de la microplaque vide doivent être dans un intervalle limité. Par exemple :  $\pm$  0,010 DO.

Si les valeurs de DO ne sont pas dans cet intervalle, il ne faut pas utiliser ce type de microplaque.

Grâce à l'utilisation de mesures de longueur d'onde doubles, l'influence de la différence dans les valeurs de DO de la microplaque est supprimée ou réduite à un niveau acceptable.



# 5.3.3 Précision de l'instrument avec des échantillons liquides

Cette procédure peut être utilisée pour vérifier la répétabilité des mesures. Il est recommandé d'utiliser une microplaque à fond plat.

Remplir une nouvelle microplaque avec une solution d'orangé G fraîchement préparée ; utiliser différentes dilutions de la solution pour chaque puits afin d'avoir un intervalle de densités optiques. Vérifier que les puits contiennent au moins 200 µl. Les séries de dilutions doivent être comprises dans l'intervalle de 0,1 à 3,0 DO. Pour parvenir à environ 3 DO, il est recommandé d'utiliser 125 mg.l<sup>-1</sup> d'orangé G (Sigma, cat. n°O7252).

Programmer un test pour utiliser le filtre à 492 nm puis mesurer la microplaque au moins trois fois.

Pour chaque puits, calculer les valeurs suivantes :

- la valeur moyenne de DO
- déviation standard

#### Exemple

#### Lectures 0,000 à 2,000 DO

La déviation standard de chaque puits doit être de (0.5 % + 0,005 OD). Calcul de la déviation maximum permise en prenant 1,000 OD comme valeur moyenne de DO :

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

#### Lectures 2,001 à 3,000 DO

La déviation standard de chaque puits doit être de (1.0 % + 0,005 OD).

Calcul de la déviation maximum permise en prenant 2,400 OD comme valeur moyenne de DO :

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

#### Lectures au-dessus de 3,000 DO

Les lectures au-dessus de 3,000 DO sont utilisées uniquement à titre indicatif et la précision ne peut être garantie.



## 5.3.4 Linéarité de l'instrument avec des échantillons liquides

Pour vérifier la linéarité de l'instrument et l'application à la longueur d'onde utilisée, utiliser la dilution sérielle d'une solution.

Le résultat dépend de la pureté du colorant utilisé et du ménisque du liquide dans les puits.

Pour référence, une dilution sérielle de solution d'orangé G pour des mesures à 492 nm peut être utilisée.

Les séries de dilutions doivent être comprises dans l'intervalle de 0,1 à 3,0 DO. Pour parvenir à environ 3 DO, il est recommandé d'utiliser 125 mg.l<sup>-1</sup> d'orangé G (Sigma, cat. n°O7252).

Pour les autres longueurs d'onde, différentes solutions peuvent être utilisées.

Pipetter alors 200 µl de chaque dilution sur la microplaque, en utilisant au moins deux échantillons pour chaque dilution afin de réduire les erreurs causées par le pipetage.

Mesurer alors la microplaque et tracer un graphique de régression linéaire de DO par rapport à la concentration à partir de la moyenne des valeurs DO mesurées.

Déterminer le carré de la valeur résiduelle non pondérée <sup>R2</sup> de la ligne de régression.

Des carrés de valeurs résiduelles typiques pour une application standard sont égaux ou supérieurs à  $R^2 = 0.998$ .



#### Note

Les données peuvent varier en raison de l'inexactitude du pipetage.



# 6. Exemple d'application

# 6.1 Introduction

Les **fichiers exemples** Magellan fournissent des méthodes et des espaces de travail Magellan afin de servir d'introduction au logiciel et de faciliter le travail des utilisateurs avec lui. Les fichiers exemples pour un dosage quantitatif et qualitatif ELISA sont installés automatiquement lors de l'installation de Magellan.

# 6.2 Exemple étape par étape : ELISA quantitatif

Ce chapitre contient un exemple étape par étape (test quantitatif) de création d'une méthode dans Magellan. En suivant les instructions, vous apprendrez comment définir des évaluations dans Magellan à partir d'une description de kit de test.



Note Les fichiers exemples apparaissent automatiquement dans la Liste des méthodes dans Magellan. Pour Magellan Tracker, ces fichiers sont disponibles dans le chemin de données par défaut, et doivent être convertis.

# 6.2.1 Description du kit de test

Les instructions suivantes se trouvent dans la description du fabricant du kit de test pour essai ELISA quantitatif IgM (détection d'anticorps) : schéma de la plaque

	1	2	3	4	5	6	7	8	9	10	11	12
A	BLK	C3	S1									
В	NC	C4	S2									
С	NC	C4	S2									
D	C1	C5	S3									
Е	C1	C5	S3									
F	C2	C6										
G	C2	C6										
Н	C3	S1										
	•	•										

BLK = blanc, NC = contrôle négatif, C1 - C6 = calibrateurs (standards), S1 - S... = échantillons



#### Mesure et évaluation

Lecture de la plaque à une longueur d'onde de 492 nm, référence à 620 nm. Sélectionnez le lecteur / plaque au puits A1 pour le blanc.

Concentration des calibrateurs (standard) :

Calibrateur 1	5 UA/ml
Calibrateur 2	10 UA/ml
Calibrateur 3	20 UA/ml
Calibrateur 4	40 UA/ml
Calibrateur 5	80 UA/ml
Calibrateur 6	160 UA/ml

Après la correction du blanc, un graphique compare les densités optiques (DO 492 - DO 620) par rapport à la concentration. La ligne de régression qui passe par ces points est la courbe standard.

Interprétation des résultats du test

IgM < 18 UA/ml	Négatif
18 UA/mI ≤ IgM < 22 UA/mI	Intermédiaire
IgM ≥ 22 UA/mI	Positif

La concentration d'IgM calculée des deux contrôles négatifs doit être inférieure à 8 UA/ml.

#### Traitement de données

Après la mesure, le fichier de données (espace de travail) est stocké automatiquement, et un rapport contenant les paramètres de mesure, le schéma de la plaque, les valeurs de blanc, la courbe standard, les concentrations IgM, la définition des valeurs seuil, les résultats qualitatifs des échantillons et les validations est créé.

De plus, le schéma et les résultats qualitatifs sont stockés dans un fichier ASCII.



# 6.2.2 Créer une méthode

Dans la Liste des assistants, sélectionnez Créer/éditer une méthode et cliquez sur OK. Cliquez sur Suivant dans la Page d'accueil de l'assistant Créer/éditer une méthode et la boîte de dialogue Sélection de fichier apparaît. Sélectionnez Créer.

Créer/éditer une méthode			
◯ Nouveau ⊙ Duvrir	Afficher:	Fichiers à partir de cet instrum 💙	Aperçu avant impression
🗀 mth	Name 🔺	Remarques	État
	<		) >
Aide Annuler <<< Retour		FAITES VOTR	E SÉLECTION
			© 2009 Tecan

#### Paramètres de mesure

Cliquez sur **Effectuer une sélection**, et la page **Paramètres de mesure** apparaît.

Paramètres de mesure - infinite F50			×
🔶 Outils laboratoire 🔹	🔷 ▼ Plaque	1 Séle	ction
<ul> <li>Plaque</li> <li>Partie de la plaque</li> </ul>	Définition de la plaque : [GRE96ft] - Greiner 96 Flat Transparent	Détails	seectorine
P Homes		Utiliser une partie de la plaque	
Absorbance	🔹 🕨 Partie de la plaque A1:H12	2	
	Absorbance	3	
Actions ×	Congueur d'onde Mesure: 405 nm V Nom: Étiquettei V		
A Déplacer plaque	Référence: 405 nm		
G Cinétique			
Cycle cinétique			
Commentaire			
Requête de l'utilisateur     Attender (barlene)			
<ul> <li>Accendre (nonoge)</li> <li>Incubation</li> </ul>			
		203	
Nombre de plaques : 1			
Aide	СНС	IISIR LES PARAMÈTRES DE LA MESURE	
Annuler			<b></b>
© 2009 Tecan			



#### 6. Exemple d'application

Dans le bandeau **Longueur d'onde**, sélectionnez 492 nm comme Longueur d'onde de mesure et 620 nm comme Longueur d'onde de référence.

🍦 🔻 Absorb	ance			3
-Longueur d'ond	e		Étiquette	
Mesure:	492 nm	*	Nom: Étiquette1 🗸	
Référence:	620 nm	*		

Continuez l'assistant en cliquant sur **Choisir les paramètres de la mesure**, et la fenêtre **Schéma de la plaque** s'affiche.

Créer/éditer une méthe Annuler Répéter Sél	o de ectionner tous les	non utilisés	10% 1	10% Zoc	m								×
Schéma de la mé	tho 🛞 🖓												
	/éditer une m	0	1	2	3	4	5	6	7	8	9	10	11 12
	éthode	A											Identificateurs
		B											ST Z 4 PC 1 0
		C											LPC Réplications - HPC Z M BF ONuméro fixe:
		D											RF Tous
		E											Rempir Supprimer sélection
		F											
		G											AN Austria
		H											P
Aide Annuler <<<	Retour												SUIVANT



#### Définir le schéma de plaque

Définissez le schéma de la plaque en utilisant la boîte de dialogue **Attribution du puits** située dans la partie droite de l'écran.



Dans la **boîte de groupe** Identificateurs, sélectionnez **BL (Blanc)**. Dans la boîte de groupe **Groupe expérimental**, le chiffre 1 reste. Dans la boîte de groupe **Répliques**, **Tous** est sélectionné automatiquement. Cliquez sur le puits **A1**, qui est alors marqué par une bordure rouge. Cliquez sur **Remplir sélection**, et le puits est étiqueté avec le type d'identificateur sélectionné.





Choisissez maintenant les réglages suivants dans la boîte de dialogue Attribution du puits :

Dans la boîte de groupe **Identificateurs**, sélectionnez **NC** (contrôle négatif). Dans la boîte de groupe **Groupe expérimental**, le chiffre **1** reste.

**Tous** est sélectionné automatiquement dans la boîte de groupe **Répliques**. En partant du puits **B1**, cliquez et glissez la souris jusqu'au **C1**. Les puits **B1** à **C1** sont alors marqués par une bordure rouge.

Cliquez sur **Remplir sélection**, et les puits sont étiquetés avec le type d'identificateur sélectionné.

Les calibrateurs (standards) doivent être assignés aux puits **D1** à **G2**. Choisissez les réglages suivants dans la boîte de dialogue **Attribution du puits** :

Dans la boîte de groupe **Identificateurs**, sélectionnez **ST (Standard)**.

Dans la boîte de groupe Expérimental, le chiffre 1 reste.

Dans la boîte de groupe Répliques, choisissez entre Numéro fixe et Tous.



#### Numéro fixe :

Seulement actifs pour les standards et les échantillons permettant d'utiliser des ID.

Si le bouton **Numéro fixe** est actif, un nombre peut être entré dans le champ de texte correspondant. Ce nombre définit combien de répliques sont prévues pour cette méthode. Le nombre de répliques entré pour chaque ID est créé dans les puits sélectionnés. Le nombre de puits sélectionné doit donc être un multiple du nombre de répliques entré.

#### Tous :

Tous les puits sélectionnés sont définis comme répliques. Si un numéro d'ID existant est choisi pour les échantillons et les standards, les puits sélectionnés sont alors ajoutés en tant que répliques aux répliques existantes. Avec tous les autres types d'identificateurs, les puits sélectionnés sont ajoutés comme répliques aux répliques existantes.

Deux boutons fléchés définissent la direction de la réplique et la séquence de numéro d'ID (horizontale ou verticale).

Pour cet exemple, sélectionnez Numéro fixe et 2.

Dans la boîte **Numéro d'ID** et dans la boîte de groupe **Répliques**, sélectionnez les **flèches verticales**.

Sélectionnez ensuite les puits D1 à G2, et cliquez sur Remplir sélection.



Note

Sélectionnez les puits comme suit : En partant du puits D1, cliquez et glissez la souris sur les puits requis jusqu'à H1. Maintenez ensuite enfoncée la touche contrôle (Ctrl), et glissez la souris sur les puits requis, d'A2 à G2.

#### Le Schéma de la plaque apparaît comme suit :



Cliquez sur **Sélectionner tous les non utilisés** dans la barre d'outils pour sélectionner tous les puits vides de la plaque. Maintenez ensuite enfoncée la touche contrôle (Ctrl) et cliquez sur le puits **H12**, afin qu'il reste vide et non marqué.



Dans la boîte de dialogue **Attribution du puis**, sélectionnez **SM (échantillon)** dans *Identificateurs*.

Dans la boîte de groupe Groupe expérimental, le chiffre 1 reste.

Dans la boîte de groupe Répliques, choisissez Numéro fixe et 2.

Laissez 1 dans la boîte **Numéro d'ID**, et dans la boîte de groupe **Répliques**, sélectionnez les **flèches verticales**. Cliquez ensuite sur **Remplir sélection**. La procédure de **Définition du schéma** est terminée.

#### Transformations

Dans la barre de contrôle située à gauche de la fenêtre, sélectionnez l'option suivante, **Ajouter nouvelle transformation...** dans l'élément **Données transformées**, afin de définir la réduction du blanc.

Une boîte de dialogue s'affiche pour vous demander si vous souhaitez définir une réduction du blanc. Cliquez sur **Oui**. La fenêtre suivante apparaît :

Créer/éditer une méthode													
Annuler Répéter		10% 100%	Zoom										
Ś	Doonées dentrée -	Donoées différer	xe.			~					ſ	Constantes	Ontions
	fx		~~					~	Données disp	onibles (séries)	te danné 🔻	Exections &Cr	ostantes V
Schéma de la plaque												1	
Valeurs de conc, dil, réf.													
Doppées transformé	1												
Réduction du blanc	0	1 9	2 3	A	5	ß	7	8	0	10	11	12	
Ajouter nouvelle transform		U 4	5 9		0		u	0					- 11
Concentrations 🛞		BL1 ST	1 3 SM1 1	SM1 5	SM1 9	SM1 13	SM1 17	SM1 21	SM1 25	SM1 29	SM1 33	SM1 37	
Courbe standard		1/1 2	2/2 2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Évaluer les données 🛞		x-BL1 x-I	BL1 x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Définition des valeurs seul		NC1 ST	1_4 SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
Validation QC	6	1/2 1 	/2 1/2 21.4 × D1.4	1/2	1/2 1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
Traitement de donn 🛞		X-DLI X-I		X-DLT	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DL1	X-DLT	X-DLT	
Exportation des données	C	NC1 SI	1_4 SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
<ul> <li>Rapport imprime</li> <li>Traitement automatisé de d</li> </ul>		x-BL1 x-I	BL1 x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
2		ST1 1 ST	1.5 SM1.3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
Divers 🛞	D	1/2 1	/2 1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
Format numérique		x-BL1 x-I	BL1 x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Notes de méthode		ST1_1 ST	1_5 SM1_3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
	E	2/2 2	12 2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
		x-BL1 x-I	3L1 x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
	E	ST1_2 ST	1_6 SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	
	F	y BI1 y	72 172 RI1 X RI1	1/2 x-RL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x-BL1	1/2 x BL1	1/2 x-BL1	1/2 x-BL1	
		STI 2 ST		SMI 8	SM1 12	SM1 16	SM1 20	SM1 24	SM1 28	SM1 32	SMI 36	SM1 40	.ep
	G	2/2 2	2/2 2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	Aust
		x-BL1 x-I	BL1 x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	AN 4
		ST1_3 SM	1_1 SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37		EO
	H	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-	3L1 X-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1		
										_			
Aide													
Ánnuler SSS Betrue											5	UIVANT	
Comment Contracted													
											© 200	9 Tecan	

**Données différence** est sélectionné automatiquement dans la boîte **Données d'entrée**. Si vous avez confirmé la définition d'une réduction du blanc auparavant, le logiciel la nomme automatiquement **Réduction du blanc** (voir Données transformées dans la barre de contrôle).

Dans la boîte **Formule**, **x-BL1** apparaît automatiquement pour cette réduction du blanc, x renvoyant à la valeur actuelle des données d'entrée dans un puits, et BL1 étant la valeur moyenne des puits blancs du groupe expérimental 1.

Pour plus de détails et d'explications concernant la définition et l'attribution des transformations, veuillez consulter la notice d'utilisation de Magellan.

Les informations suivantes apparaissent dans chaque puits (exemple : puits A5) :

SM1_9	Echantillon, groupe expérimental numéro 1, ID d'échantillon numéro 9.
2/2	Le numéro de réplique est 2, le nombre total de répliques est 2.
x-BL1 ou 1	Transformation x-BL1 assignée (lorsque Transformation est sélectionné) ou valeur de facteur de dilution de 1 (lorsque Valeurs conc., dil., réf. est sélectionné).



#### Définition de valeur de concentration / dilution / référence

Dans la barre de contrôle, sélectionnez **Valeurs conc., dil., réf.** dans l'élément **Schéma de la méthode** pour définir les valeurs respectives en suivant la description du kit de test.

Calibrateur 1	5 UA/ml
Calibrateur 2	10 UA/ml
Calibrateur 3	20 UA/ml
Calibrateur 4	40 UA/ml
Calibrateur 5	80 UA/ml
Calibrateur 6	160 UA/ml

Assurez-vous que **ST** est sélectionné dans la liste **Sélectionner identificateur**.

Dans la liste **Identificateur**, une liste des standards du groupe exp. 1 apparaît. Dans la boîte **Concentration** correspondante de **ST1\_1**, entrez le nombre **5**, et dans la boîte **Unité**, tapez UA/ml. Dans la boîte **Concentration** correspondante de **ST1\_2**, tapez le nombre **10**. L'unité n'a besoin d'être définie qu'une seule fois, et est valide pour tous les standards. Entrez les valeurs pour ST1\_3 à ST1\_6 de la même manière.

L'écran montrant le schéma de la plaque et la concentration s'affiche :





#### **Courbe standard**

Dans la barre de contrôle, cliquez sur **Courbe standard** dans l'élément **Concentrations** pour définir la courbe standard appropriée.

On trouve ce qui suit dans la description du kit de test :

Après la correction du blanc, un graphique compare les densités optiques (DO 492 - DO 620) par rapport à la concentration. La ligne de régression qui passe par ces points est la courbe standard.

Dans l'onglet **Données**, sélectionnez **Réduction du blanc** comme données d'entrée.

Données Type d'ana	lyse Interception Axe Graphique
Données d'entrée :	Réduction du blanc
<ul> <li>Standards</li> </ul>	du schéma
O Standards fichier exté	venant du Sélectionner
O Standards expériment	venant du groupe
🔿 Pas de trac	té de la courbe standard
Concentration	ns supplémentaires

Dans l'onglet Type d'analyse, sélectionnez Régression linéaire.

<ul> <li><u>Régression linéaire</u></li> <li><u>Régression non linéaire</u></li> </ul>		Echelle des données :	
○ Spline cubique ○ Akima		endolend) A	
O Polynôme	2		
O Quatre Paramètres			
🔘 Quatre Paramètres Marquardt		Dha	
🔘 Cinq Paramètres		Plus	



### 6. Exemple d'application

Dans l'onglet **Axe**, définissez la légende et l'échelle des axes comme cidessous :

Libelle : Concentration [U	A/mlj		
Couleur :		Echelle logarithmique	
<ul> <li>Sélectionner automatiquem</li> </ul>	ent l'étendue		
○ Étendue	min. :	max,:	
🔽 Quadrillage	Couleur :	▼ Style :	
ixe Y			
Libellé : Réduction du blan	ic		
Couleur :		Echelle logarithmique	
<ul> <li>Sélectionner automatiquem</li> </ul>	ent l'étendue		
OEtendue	min. :	max.:	
🗹 Quadrillage	Couleur :	🖌 Style :	

Dans l'onglet **Graphique**, définissez le titre du graphique, les courbes, la police et l'affichage du graphique.

Libellé :	IgM-ELISA			
Couleur :	•			
Courbes				
		Libellé: Grp. 1		
Couleur	•	Masquer la courbe		
Symbole :	<b>—</b>	Largeur de ligne 🛛 🗧		
Taille :	moyen 💌			
Police		Afficher		
💿 petit		🔽 Légende	Interception	
O moyen		Points de base	Barres d'erreurs	



#### Définir les valeurs seuil

Dans la barre de contrôle, sélectionnez **Définition des valeurs seuil** dans l'élément **Evaluer les données** pour définir les limites de l'évaluation qualitative.

La description du kit de test contient les instructions suivantes :

Interprétation des résultats du test :

IgM < 18 UA/mI	Négatif
18 UA/ml ≤ IgM < 22 UA/ml	Intermédiaire
IgM ≥ 22 UA/mI	Positif

Utilisez la procédure suivante pour définir les valeurs seuil appropriées :

Dans la boîte Données d'entrée, sélectionnez Conc. moyenne (UA/ml).

Le tableau **Valeurs seuil** représente une échelle indiquant les bornes supérieures et inférieures des **Limites** et des **Etiquettes**. Dans **Limites**, tapez 22 comme première limite (supérieure) et 18 comme seconde limite (inférieure).

Dans **Etiquettes**, entrez l'interprétation du test (**Positif**, **Intermédiaire**, et **Négatif**) dans les boîtes individuelles. Utilisez la palette de couleur déroulante pour assigner une couleur :

- Positif rouge
- Intermédiaire bleu
- Négatif vert

Le contenu de l'écran est le suivant :

Données d'entrée :	Conc.moyenne. (UA/ml)	
Valeurs seuil Couleurs El	Cont. Hoyenne. (CA)/III)       Itiquettes       Itiquettes	
Saisie de formule Variable BL1	Opérateurs Fonctions	Test compétitif

Cliquez sur **Sélection des résultats de valeurs seuil** pour sélectionner les types d'identificateurs pour lesquels les résultats de valeurs seuil doivent être affichés.

#### Définir les validations QC

Dans la barre de contrôle, cliquez sur **Validations QC** dans l'élément **Evaluer les données**. Des critères de validation doivent être définis pour le test, de manière à garantir la validité des résultats du test.

Dans cet exemple, la condition suivante doit être remplie :

La concentration d'IgM calculée des deux contrôles négatifs doit être inférieure à 8 UA/ml.

Dans la boîte Données d'entrée, sélectionnez **Conc. unique (UA/ml)**. Dans la première ligne, tapez **NC1\_1<8**.



Note NC1\_1 signifie Contrôle négatif du groupe expérimental 1, réplique 1.

Dans la seconde ligne, tapez NC1\_2<8.

La boîte de dialogue Validations QC s'affiche alors ainsi :

Conditions do vol	dation			
NC1 1-9	dation			
NC1_1<8				
1				
Saisie de formule				
Variable	Onérateur	s Fonctions		
C Plaque à plaque	✓ … +	Y and	×	
Données d'entrée : [	+			
BL1 CC Plaque à plaque Données d'entrée : [ Contrôle 1:	Données différence	Moyenne :	s:	Ť
Données d'entrée : [ Contrôle 1: Contrôle 2:	Données différence	Moyenne :	s:	Ť
Données d'entrée : [ Contrôle 1: Contrôle 2: Contrôle 3:	Données différence	Moyenne : Moyenne : Moyenne	s:           s:           s:	Ť
BL1 CC Plaque à plaque Données d'entrée : [ Contrôle 1: Contrôle 2: Contrôle 3: Contrôle 4:	Données différence       BL1       BL1       BL1       BL1	Moyenne : Moyenne Moyenne Moyenne :	S:           S:           S:           S:           S:           S:	↑ ↓
Données d'entrée : [ Contrôle 1: Contrôle 2: Contrôle 3: Contrôle 4:	<ul> <li>Connées différence</li> <li>BL1 </li> <li>BL1 </li> <li>BL1 </li> <li>BL1 </li> <li>SL1 </li> </ul>	Moyenne : Moyenne Moyenne Moyenne :	s:       s:       s:       s:	↑ ↓
Données d'entrée : [ Contrôle 1: Contrôle 2: Contrôle 3: Contrôle 4: Otrous les espace Espaces de trav		Moyenne : Moyenne Moyenne Moyenne : miers 1 jours	s:       s:       s:       s:	↑ ↓



# 6.2.3 Organiser le rapport imprimé

Dans la barre de contrôle, cliquez sur **Rapport imprimé** dans l'élément **Traitement de données**. L'écran suivant s'affiche :

⊙ Liste ○ Matrice	Vous pouvez utiliser le glisser-déposer pour créer un rapport. Choisir un élément dans la liste des données disponibles et le faire glisser dans la liste des données sélectionnées. Le déposer à la position appropriée dans le rapport.
onnées disponibles:	Données sélectionnées:
	Insérer Ajouter  E Collection données Haut Bas Haut Bas Propriétés Supprimer to
E Signature	Tmoster

Dans l'onglet **Sélection de données**, toutes les données de rapport disponibles sont contenues dans la boîte **Données disponibles**. En utilisant les boutons **Insérer et Ajouter**, les données peuvent être transférées dans la boîte **Données sélectionnées**. Les données peuvent également être transférées par glisserdéposer.

Dans la boîte **Imprimer en tant que**, choisissez d'imprimer les données sous forme de matrice ou sous forme de liste avec une orientation spécifique.

Dans cet exemple, un rapport contenant les paramètres de mesure, le schéma de la plaque, les valeurs de blanc, la courbe standard, les concentrations IgM, la définition des valeurs seuil, les résultats qualitatifs des échantillons et des validations devrait être créé.

Avant de créer le rapport, l'option par défaut Liste verticale/Données différence doit être retirée de la boîte Données sélectionnées. Seul Paramètres de mesure doit rester dans la boîte Données sélectionnées. Imprimer en tant que liste doit être changé en Imprimer en tant que matrice.



#### 6. Exemple d'application

Liste   Matrice	Vous pouvez ut dans la liste de sélection	iliser le glisser-déposer pour créer un rapport s données disponibles et le faire glisser dans nées. Le déposer à la position appropriée da	. Choisir un élément la liste des données ns le rapport.
onnées disponibles:		Données sélectionnées:	
<ul> <li>Données différence - Déviati ▲</li> <li>Données transformées</li> <li>Concentrations</li> <li>Concentrations</li> <li>A Résultats qualitatifs</li> <li>'s, ID d'échantillon</li> <li>Schéma de la méthode</li> <li>(rithers de validation QC</li> <li>Paramètres de mesure</li> <li>R Remarques</li> <li>Protocole d'erreur</li> <li>Trace d'audit</li> <li>Signature</li> <li><saut de="" page=""></saut></li> <li><séparateur></séparateur></li> <li><séparateur></séparateur></li> <li><séparateur></séparateur></li> <li><séparateur></séparateur></li> <li><séparateur></séparateur></li> <li><séparateur></séparateur></li> </ul>	Insérer Ajouter	■ E. Collection données           □	Haut Bas Propriétés Supprimer Supprimer tout
	Importer		

Sélectionnez Schéma de méthode/Schéma dans la boîte Données disponible, et attachez-le au rapport en cliquant sur Ajouter. Ensuite, insérez Réduction du blanc, Conc. moyenne (UA/ml) et Résultats de valeur seuil dans la matrice en sélectionnant les éléments correspondants et en cliquant sur Insérer.

Ajoutez Graphique : Courbe standard, définition de valeurs seuil et Critères de validation QC aux données sélectionnées. L'étape de paramétrage des données de la procédure de définition du rapport est terminée ; la boîte de dialogue Rapport imprimé ressemble à ceci :



Dans les onglets **En-tête** et **Pied de page**, définissez la mise en page de l'entête et du pied de page du rapport (veuillez consulter la notice d'utilisation de Magellan pour plus de détails).





#### Exportation de données

Dans la barre de contrôle, cliquez sur **Exportation de données** dans l'élément **Traitement de données**. Dans cet exemple, le schéma et les résultats de valeurs seuil doivent être enregistrés dans un fichier ASCII. Sélectionnez le **Schéma** et les résultats de **Valeurs seuil** dans la fenêtre **Données disponibles** ; cliquez sur la flèche → pour les insérer dans la fenêtre **Données sélectionnées**. L'écran affiche alors les informations suivantes :

Données disponibles :			Données sélectionnées :	
🖅 🛍 Données instrument	^		Schéma	Haut
🖨 🤆 Donnée réduite			Donnees difference	- Read
Données différence				Das
Données différence - Moyen				
Données différence - Déviation sta				
Données différence - Coefficient de				
🕀 🔏 Données transformées				
😥 👔 Concentrations		->		
Résultats de valeur seuil				
		<u></u>		
🖃 🛲 Schéma de la méthode				
Positions de puits				
Noms de la méthode sur bande	-			
Concentrations originales				
Facteurs de dilution	~			
	-			
		1	·]	
Options d'exportation	r AS	5CII)	Exporter dans Excel	
() ()				



Note Les données exportées doivent toujours contenir le schéma ou la liste ID d'échantillons.

#### Traitement automatisé de données

Dans la barre de contrôle, cliquez sur **Traitement automatisé de données** dans l'élément **Traitement de données**.

Automatique	
Chargement liste ID d'échantillon	Plus
Enregistrer l'espace de travail	Plus
Exporter dans un fichier ASCII	
Exporter dans un fichier ASTM (LIS)	Plus
Exporter dans Excel	
Imprimer	
Afficher les résultats après mesures	Plus

Sélectionnez **Exporter dans un fichier ASCII**, et **Afficher les résultats après mesures**. Dans Magellan **Tracker**, **Enregistrer l'espace de travail** est sélectionné par défaut, et ne peut pas être modifié.



### Enregistrer la méthode

Cliquez sur **Suivant** pour ouvrir la fenêtre **Enregistrer sous**. Entrez le nom de fichier de la méthode et remplissez tout autre champ si nécessaire.

Créer/éditer une	e méthode			
Enregistrer sous:	mth	Name 🔺	Remarque:	État
		Ivanic —	rremarques	
		<		>
Nom du fichier:	Méthode1.mth			
Remargues fichier:				
				~
Commenter trace d'a	audit			
				~
Organiser les	favoris Trace d'audit Si	gnatures Mot de	passe de la méthode:	
Aide	Dér	narrer cette méthode maintenant	ENDECISTDED	
Annuler	<	and settle more and manyoridin	TERMINEI	Ř 🔽
			© 2009 Tecar	1

Champ de texte Nom de fichier	Un nom de fichier doit être entré. Un nom de fichier par défaut est suggéré automatiquement, mais il peut être modifié.
Champ de texte Remarques fichier	Les commentaires entrés ici seront enregistrés et affichés avec le nom du fichier.
Champ de texte Commenter trace d'audit	Les commentaires entrés ici seront enregistrés dans la trace d'audit. Cette option n'est disponible qu'avec <b>Magellan Tracker</b> .
Bouton Organiser les favoris	La boîte de dialogue <b>Organiser les favoris</b> apparaît.
Champ de texte Mot de passe de la méthode	Entrez un mot de passe pour protéger la méthode.
Case à cocher Démarrer cette méthode maintenant	La méthode sera démarrée immédiatement après avoir cliqué sur <b>Enregistrer&amp;Terminer</b> .

Pour plus de détails, veuillez consulter la notice d'utilisation de Magellan.



# 6.2.4 Exécuter la méthode

Si **Démarrer cette méthode maintenant** est sélectionné dans la boîte de dialogue **Enregistrer sous** de **l'Assistant Créer/éditer une méthode**, la boîte de dialogue **Assistant Démarrer la mesure/Démarrer la mesure** apparaîtra après avoir cliqué sur **Enregistrer**. Dans la boîte de dialogue **Démarrer la mesure**, un nom d'espace de travail par défaut est assigné. Ce nom peut être modifié par l'utilisateur s'il le souhaite.

Démarrer mesure	
Mesure         Espace de travail:       05112009-001.wsp         Méthode:       Méthode1.mth         Liste des ID d'échantillon:       Insérer         Instrument       Utiliser dispositif d'empilement       Actuelle:         Utiliser dispositif d'empilement       Cible:       n. déf         Plaque rentrée       Contrôle de temp       Cible:       n. déf         Mouvements       Contrôle injecteur       Optimise la position Z	Paramètres de mesure Plaque Description de plaque: [GRE96it] - Greiner S Plaque avec couvercle: Non Code barre: Non Partie de la plaque Etendue: A1:H12 Absorbance Longueur d'onde de mesure: 492 nm Longueur d'onde de référence: 620 nm Étiquette: Étiquette1
Aide Annuler <<< Retour	DÉMARRER
	© 2009 Tecan

Cliquez sur **Démarrer** pour lancer la mesure. Un espace de travail sera créé automatiquement, contenant toutes les informations entrées auparavant, et qui collectera toutes les valeurs de mesure. Pendant que la mesure est exécutée, une boîte de dialogue de statut de la mesure apparaît, pour indiquer la progression de la mesure.

Lorsque la mesure est terminée, la boîte de dialogue **Résultats** apparaît, dans laquelle tous les résultats et calculs peuvent être consultés. Des messages d'erreurs peuvent s'afficher lorsqu'une mesure est effectuée sans les liquides correspondants (par ex. standards).



## 6.2.5 Évaluer le résultat

Sélectionnez **Évaluer les résultats** pour voir et évaluer les données brutes. Les paramètres d'évaluation peuvent être visualisés et les données réévaluées.

Cette section vous guide tout au long de l'**Assistant Évaluer les résultats**, en utilisant un exemple de fichier d'espace de travail installé automatiquement lors de l'installation de Magellan.

Dans la boîte de dialogue Liste des assistants, cliquez sur Évaluer les résultats.

Cliquez sur **Suivant** dans la **Page d'accueil** de **l'assistant Évaluer les résultats** et la boîte de dialogue **Sélection de fichier** apparaît.

Sélectionnez l'espace de travail Quantitative ELISA

**example\_Sunrise\_InfiniteF50.wsp** dans la liste de fichiers et cliquez sur **Faites votre sélection**. Les calculs sont exécutés et la fenêtre de schéma de la plaque est affichée :

Évaluer les résultats - Quantit	ative l	ELISA example	_Sunrise_Infi	initeF50.ws	o - Quantitat	live ELISA e	cample_Sun	rise.mth							×
Fichier Éditer Instrument															
Editer			10% 100%	Zoom											
	0														
lu.															$\sim$
Données instrument 🛞	Éval	0	4	2	2	A	B	ß	7	8	0	10	11	12	
Données de référence	ler		U	6	0		0	U	U	0	0	10	0.0	025	1
Données de mesure - Moyen	C S D		DI 1	CT1 2	Ch41_4	CMI 5	Ch41 0	CM1 12	CM41 47	CL41 . 24	CM1 25	CM41 20	CM1 22	CM1 27	
s - Données de mesure	Ésul		1/1	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
v - Données de mesure	ats	6-3	0 004	0 207	0.1	0.816	0 174	0 166	0.083	0.085	0.085	0.08	0.162	0.131	
Données de référence - Mo			NICI	OT4 4	CL41_2	Ch41_C	Ch41 40	CM4 44	CM4 40	CM4 00	CM4	CM4 20	CM4 . 24	Ch41 20	
s - Données de reference y - Données de référence		R	1/2	311_4	SIVIT_2	SIVIT_0	SIVIT_10	3WI 14	3NII_10	3WI _22	SIVIT_20	5IVIT_50	5WI1_04	31/2	
Données de mesure comme	Édit		0.060	0.418	0.784	0.212	0.106	0.156	0.106	0.216	0.123	0 103	0.131	0.152	
Données de référence com	- P		0.000	0.410	0.704	0.212	0.100	0.100	0.100	0.210	0.120	0.100	0.101	0.102	
	tho	R	2/2	511_4	SIVIT_2	SIVI1_6	301_10	SIVI1_14	31/11_18	3MI_22	31/1_26	301_30	3MT_34	3101_38	
Donnee reduite	6		2/2	2/2	0.764	0.205	2/2	0.155	0.104	0.24	0.120	0.112	0.125	212	
Données différence - Moyen			0.000	0.410	0.704	0.203	0.195	0.133	0.104	0.21	0.125	0.112	0.123	0.133	
s - Données différence		D	511_1	511_5	SM1_3	SM1_/	SM1_11	SM1_15	SM1_19	SM1_23	SM1_2/	SM1_31	SM1_35	SM1_39	
v - Données différence			0.052	0.000	1/2	0.000	1/2	0.400	0.424	0.44	0.460	0.405	0.427	1/2	
Données différence comme			0.052	0.030	0.04	0.065	0.096	0.106	0.121	0.11	0.102	0.105	0.127	0.095	
Données transformé(♡)		E	ST1_1	SI1_5	SM1_3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
2		E	212	212	2/2	212	2/2	212	2/2	212	2/2	212	2/2	2/2	
Concentrations 🛞			0.051	0.84	0.629	0.085	0.1	0.11	0.125	0.115	0.100	0.111	0.129	0.09	
Résultats qualitatifs (😒		l e	S11_2	SI1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	
11 A		F	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
Schéma de la métho 📎			0.103	1.000	0.323	0.104	0.076	0.155	0.145	0.105	0.112	0.094	0.155	0.145	10
Validation QC 🛞		B	ST1_2	ST1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	Istri
Critères de validation QC		G	2/2	ZIZ	212	2/2	2/2	2/2	2/2	2/2	212	2/2	2/2	2/2	4 At
			0.103	1.000	0.314	0.099	0.079	0.155	0.130	0.104	0.110	0.092	0.124	0.149	CAP
Remarques		1 10	ST1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37		μ
Protocole d'erreur		l li	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2		JI
Trace d'audit			0.206	0.105	0.845	0.199	0.167	0.085	0.081	0.082	0.081	0.167	0.127		
Signature														_	
		Paramètres de r	mesure												1
		SUNRISE													
		Description de p	laque: [SUNRISE,	_Standard] - Gr	einer 96 Flat										3
Aide															
	1												:	SUIVANT	
Aunuer <<< Retour	1														
													@ 200	O Tarres	

La valeur calculée est affichée dans chaque puits individuel. La fenêtre de schéma de la plaque change en fonction de l'élément sélectionné dans la barre de contrôle. Les paramètres et les réglages peuvent être modifiés en utilisant les éléments de la barre de contrôle. Si la méthode doit être modifiée, cliquez sur l'onglet **Editer méthode**.



Cliquez sur le puits avec le bouton droit de la souris pour faire apparaître le menu contextuel :

SM	1_10	SM1	_14	SM1	_18	SN	
2 0.1	<b>Résu</b> Déta Édite	Résumé Détails Éditer					
SM 1 0.(	Édite Copi Colle	Éditer paramètres cinétiques Copier les paramètres cinétiques Coller les paramètres cinétiques					
SM 2 0	Graphique: multi-étiquetage Graphique: cinétique Graphique: multi-étiquetage Graphique: courbes FLT						
SM	Grap Grap	Graphique: Spectra Graphique: Séries de dilutions					
0.0	Masquer/démasquer la sélection Afficher/Masquer la mise en page						

En sélectionnant **Résumé**, la fenêtre suivante s'affiche et fournit des informations détaillées sur la définition et les réglages du puits sélectionné :

Puits: C5					
📮 👜 Schéma de la méthode					
🖅 🛲 Identificateur					
🗉 👜 Alias de liquide					
Đ 📠 Repliquer					
🕀 📠 Dilution					
🗊 🛲 Réduction données multi-étiquetage					
🛨 📠 Formule(s) de transformation					
Formule(s) de transformation cinétique					
Formule(s) de transformation de concentration	- Eormule(s) de transformation de concentration				
🕀 👜 Noms de la méthode sur bande					
🕣 🛄 Données instrument					
🗄 💥 Donnée réduite					
🗉 🔟 Moyenne					
🗉 🔏 Données transformées					
Concentrations					
🗄 🔟 Statistiques					
Gauche Haut Droite Bas					
vévelopper tou	ОК				

Cliquez sur **Suivant** dans la fenêtre de schéma de la plaque et la boîte de dialogue **Enregistrer sous** apparaît. Vous pouvez y entrer un nom de fichier et des remarques. Cliquez sur le bouton **Enregistrer** situé à gauche de la fenêtre pour enregistrer le fichier ; vous pouvez continuer à travailler avec la méthode ou l'espace de travail. Cliquez sur le bouton **Terminer** situé en bas à droite de l'écran pour enregistrer le fichier et fermer l'assistant. Le programme retourne alors à la liste des assistants.



## 6.2.6 Résumé de la définition de l'ELISA quantitatif dans Magellan

#### 1. Soustraire la valeur du blanc

#### Définitions dans Magellan

Cliquez sur **Ajouter nouvelle transformation** dans la barre de contrôle ; une fenêtre apparaît où l'on vous demande si vous voulez définir une **Réduction du blanc**. Cliquez sur **Oui** et la formule de **Réduction du blanc** est assignée automatiquement à tous les puits.

#### 2. Définir les concentrations

# Définitions dans Magellan (barre de contrôle – Schéma de méthode / Valeurs conc., dil., réf.)

Sélectionnez l'identificateur : ST

Unité : UA/ml

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

#### 3. Définir la courbe standard

# Définitions dans Magellan (barre de contrôle – Concentrations / Courbe standard)

Données d'entrée réduction du blanc

Type d'analyse	régression linéaire
Axe X	linéaire
Axe Y	linéaire

#### 4. Définir les valeurs seuil

#### Définitions dans Magellan (barre de contrôle – Évaluer les données / Définir les valeurs seuil)

Données d'entrée :Conc. moyenne (UA/ml) Limites : 22 18 Positif ≥ 22 > intermédiaire ≥ 18 > négatif Test non compétitif

#### 5. Validation QC

# Définitions dans Magellan (barre de contrôle – Évaluer les données / Validation QC)

Données d'entrée : Conc. unique (UA/mI)

Condition de validation 1NC1\_1<8

Condition de validation 2NC1\_2<8

NC1\_1.....Contrôle négatif première réplique premier groupe expérimental

NC1\_1.....Contrôle négatif deuxième réplique premier groupe expérimental



# 7. Nettoyage, maintenance et élimination des déchets

# 7.1 Introduction

Ce chapitre contient les procédures suivantes :

- Nettoyer l'instrument
- Désinfecter l'instrument
- Assurer la maintenance de l'instrument
- Insérer ou remplacer des filtres dans la roue porte-filtres
- Instructions d'élimination





AVANT TOUTE OPÉRATION DE NETTOYAGE OU DE MAINTENANCE, RETIREZ LA MICROPLAQUE.



### AVERTISSEMENT

DÉBRANCHEZ L'INSTRUMENT DE L'ALIMENTATION EXTERNE AVANT LE NETTOYAGE ET LA DÉSINFECTION.



#### PRÉCAUTION

NE DÉPLACEZ PAS LE TRANSPORT DE PLAQUE MANUELLEMENT, SAUF SI L'INSTRUMENT EST HORS TENSION.

# 7.2 Nettoyer l'instrument



### AVERTISSEMENT

LA PROCÉDURE DE NETTOYAGE DOIT ÊTRE EFFECTUÉE DANS UNE PIÈCE BIEN VENTILÉE PAR DU PERSONNEL FORMÉ ET AUTORISÉ, PORTANT DES GANTS À USAGE UNIQUE ET DES LUNETTES ET DES VÊTEMENTS DE PROTECTION.

Nettoyez le boîtier de l'instrument et le transport de plaque uniquement avec un chiffon sec ou humide. S'il est très sale, nettoyez l'instrument avec un chiffon imprégné d'une solution à 70% maximum d'éthanol ou un détergent doux, du Microcide SQ, ou du Decon 90. Essuyez avec une serviette en papier.

Si un liquide quelconque est renversé sur l'instrument, il doit être essuyé immédiatement pour éviter qu'il n'atteigne le système optique et ne cause une perte de performances ou une erreur.



# 7.3 Désinfection de l'instrument



#### AVERTISSEMENT

#### SI UN LIQUIDE RENVERSÉ SUR LE TRANSPORT DE PLAQUE EST POTENTIELLEMENT INFECTIEUX, IL DOIT ÊTRE DÉSINFECTÉ SELON LES LOIS ET RÉGULATIONS NATIONALES APPLICABLES.

Toutes les parties de l'instrument qui sont entrées en contact avec des échantillons biologiques, des échantillons de patients, des échantillons de contrôle positifs ou des matériels dangereux doivent être traitées comme des zones potentiellement infectieuses.



#### AVERTISSEMENT LA PROCÉDURE DE DÉSINFECTION ET LES DÉSINFECTANTS DOIVENT ÊTRE EN CONFORMITÉ AVEC LES LOIS ET LES RÉGULATIONS NATIONALES APPLICABLES.



#### AVERTISSEMENT

IL EST TRÈS IMPORTANT QUE L'APPAREIL SOIT SOIGNEUSEMENT DÉSINFECTÉ AVANT DE LE SORTIR DU LABORATOIRE OU AVANT QU'UNE OPÉRATION DE MAINTENANCE SOIT ENTREPRISE.

Avant de rendre l'instrument au distributeur local ou à un centre d'assistance, il faut désinfecter toutes les surfaces et le transport de plaque, et l'autorité opérationnelle doit remplir un certificat de sécurité. Si le certificat de sécurité n'est pas fourni, le distributeur local ou le centre d'assistance peuvent refuser l'instrument, ou les douanes peuvent décider de le retenir.

## 7.3.1 Solutions de désinfection

Les surfaces externes et le transport de plaque de l'instrument doivent être désinfectés à l'aide d'une solution désinfectante telle que :

- Microcide SQ
- Decon 90
- Ethanol à 70%



RISQUE D'INCENDIE ET D'EXPLOSION ! LES ALCOOLS TELS QUE L'ÉTHANOL OU L'ISOPROPANOL SONT INFLAMMABLES ET PEUVENT CONDUIRE À DES EXPLOSIONS ET/OU DES INCENDIES LORSQU'ILS SONT TRAITÉS DE FAÇON INCORRECTE. DES PRÉCAUTIONS DE SÉCURITÉ DE LABORATOIRE ADAPTÉES DOIVENT ÊTRE RESPECTÉES.

**AVERTISSEMENT** 



PRÉCAUTION N'UTILISEZ JAMAIS D'ACÉTONE, CELA ENDOMMAGERAIT LES COUVERCLES.



## 7.3.2 Procédure de désinfection

Si le laboratoire ne possède pas de procédure de désinfection spécifique, suivez la procédure suivante pour désinfecter les surfaces externes et le transport de plaque de l'instrument.







## AVERTISSEMENT

LA PROCÉDURE DE DÉSINFECTION DOIT ÊTRE EFFECTUÉE DANS UNE PIÈCE BIEN VENTILÉE PAR DU PERSONNEL FORMÉ ET AUTORISÉ, PORTANT DES GANTS À USAGE UNIQUE ET DES LUNETTES ET DES VÊTEMENTS DE PROTECTION.

#### PRÉCAUTION

LE DÉSINFECTANT DE SURFACE PEUT EXERCER UNE INFLUENCE NÉGATIVE SUR LES PERFORMANCES DE VOTRE INSTRUMENT S'IL EST UTILISÉ À L'INTÉRIEUR DE CELUI-CI OU S'IL Y PÉNÈTRE ACCIDENTELLEMENT.

#### AVERTISSEMENT

DÉBRANCHEZ L'INSTRUMENT DE L'ALIMENTATION PRINCIPALE AVANT DE LE DÉSINFECTER POUR ÉVITER TOUT RISQUE D'INCENDIE OU D'EXPLOSION.

- 1. Portez des gants, des vêtements et des lunettes de protection.
- 2. Préparez un conteneur approprié pour tous les matériels jetables utilisés pendant la procédure de désinfection.
- 3. Déconnectez l'instrument du logiciel et mettez-le hors tension.
- 4. Sortez soigneusement le transport de plaque de l'instrument.



- Appliquez soigneusement le désinfectant sur le transport de plaque en suivant les instructions du fabricant. N'utilisez pas trop de désinfectant pour éviter que la solution ne pénètre dans l'instrument ou ne se dépose sur les lentilles lorsque vous rentrerez le transport de plaque dans l'appareil.
- Après avoir laissé agir le produit (pendant la durée préconisée par le fabricant dans ses instructions), essuyez le transport de plaque avec une serviette en papier douce et un détergent doux ou de l'eau distillée pour enlever toute trace de désinfectant.
- 7. Rentrez soigneusement le transport de plaque de l'instrument.



7. Nettoyage, maintenance et élimination des déchets



- 8. Appliquez soigneusement le désinfectant sur la plaque de base du transport de plaque.
- 9. Après avoir laissé agir le produit, essuyez la plaque de base du transport de plaque avec une serviette en papier douce et un détergent doux ou de l'eau distillée pour enlever toute trace de désinfectant.
- 10. Appliquez soigneusement la solution désinfectante sur toutes les surfaces externes de l'instrument.
- 11. Après avoir laissé agir le produit, essuyez l'instrument avec une serviette en papier douce et un détergent doux ou de l'eau distillée pour enlever toute trace de désinfectant.
- 12.Essuyez les surfaces externes de l'instrument avec une serviette en papier douce.
- 13.Renouvelez la procédure de désinfection sur tous les accessoires déplacés ou renvoyés.
- 14. Eliminez le conteneur avec le matériel jetable selon les lois et les régulations nationales applicables.
- 15. Désinfectez vos mains et nettoyez-les avec un détergent doux.
- Lorsque vous renvoyez l'instrument au distributeur local/centre d'assistance, veuillez suivre les étapes suivantes :
- 16.Emballez l'instrument et ses accessoires.
- 17.Remplissez le certificat de sécurité (voir ci-dessous) et fixez-le sur l'extérieur de la boîte, de façon visible.

## 7.3.3 Certificat de sécurité

Afin d'assurer la sécurité et la santé du personnel, il est demandé à nos clients de remplir un **Certificat de sécurité** (livrée avec l'instrument) et d'en fixer une copie sur le dessus de l'emballage dans lequel l'instrument est renvoyé (visible de l'extérieur de l'emballage d'expédition !), et de joindre une seconde copie aux documents d'expédition avant de l'envoyer au centre d'assistance pour révision ou réparation.

L'instrument doit être désinfecté sur le site de l'autorité opérationnelle avant son expédition (voir 7.3.2 Procédure de désinfection).

La procédure de désinfection doit être effectuée dans une pièce bien ventilée par du personnel autorisé et formé, et portant des gants non poudrés à usage unique et des lunettes et des vêtements de protection.

La procédure de désinfection doit être effectuée conformément aux régulations nationales, régionales et locales.

Si le certificat de sécurité n'est pas fourni, le centre d'assistance peut refuser l'instrument

Votre représentant Tecan local peut vous faire parvenir une nouvelle copie du certificat de sécurité, si nécessaire.



# 7.4 Plan de maintenance préventive pour INFINITE F50 PLUS

Les procédures de maintenance préventive suivantes sont recommandées.

## 7.4.1 Tous les mois

Nettoyer le boîtier et le transport de plaque avec un détergent doux au moins une fois par mois, plus souvent si nécessaire.



#### PRÉCAUTION N'UTILISEZ JAMAIS D'ACÉTONE, CELA ENDOMMAGERAIT LES COUVERCLES.

# 7.4.2 Tous les quatre ans

Il est recommandé de remplacer les filtres tous les 4 ans.

# 7.5 Remplacement et installation du filtre

L'INFINITE F50 PLUS doit être connecté à Magellan pour effectuer une insertion ou un remplacement de filtre guidé par le logiciel. Si la connexion est perdue au cours de la procédure, en raison d'une interruption accidentelle de la connexion entre l'instrument et l'ordinateur, Magellan doit être arrêté et l'instrument doit être mis hors tension. Si cela se produit, continuez la procédure comme indiqué cidessous. Lorsque vous aurez terminé, rétablissez la connexion en redémarrant l'appareil et Magellan, et définissez les filtres nouvellement insérés.



#### PRÉCAUTION LORS DE LA MANIPULATION DES FILTRES, FAITES ATTENTION À NE PAS LES ENDOMMAGER OU LES SALIR AVEC DES EMPREINTES DE DOIGT OU DE LA POUSSIÈRE.



## 7.5.1 Procédure de changement de filtre

Les filtres de la roue porte-filtres standard peuvent être remplacés ou complétés en suivant la procédure suivante :

- 1. Dans la fenêtre de liste des assistants, cliquez sur Divers.
- 2. Cliquez sur Contrôle instrument
- 3. Cliquez sur Définir porte-filtres...
- 4. Cliquez sur Changement de filtre pour lancer la procédure.
- 5. Retirez toute microplaque du transport de plaque !
- 6. Penchez soigneusement l'instrument vers l'arrière jusqu'à ce qu'il repose sur son côté arrière, le fond de l'instrument face à vous.
- 7. Retirez la plaque de protection du dessous de l'appareil en dévissant les vis orange.



8. Retirez la roue porte-filtres à fixation magnétique en la sortant soigneusement de l'instrument.



9. Placez la roue porte-filtres sur une surface plane et propre.



10.Pour remplacer un filtre, utilisez l'outil de montage de filtre pour retirer le filtre de son logement.

Veuillez contacter votre distributeur local Tecan pour plus d'informations sur l'outil de montage de filtre et les filtres disponibles.



- 11. Alignez l'outil de montage de filtre avec l'encoche de la bague de rétention. Tournez l'outil et retirez la bague de rétention en la tirant hors du logement du filtre.
- 12. Retournez la roue porte-filtres de manière à faire glisser le porte-filtres hors du logement. N'utilisez pas l'outil de montage pour pousser les filtres hors de leur logement, car cela pourrait endommager le filtre.
- 13.Un nouveau filtre doit être inséré dans le logement dans le bon sens, et en veillant à ne pas endommager le filtre, ni à mettre d'empreintes de doigt dessus.





Note Assurez-vous que le filtre est inséré correctement.



#### 7. Nettoyage, maintenance et élimination des déchets

14.Placez la bague de rétention à l'extrémité de l'outil de montage de filtre et tournez-la de manière à ce qu'elle ne puisse pas glisser.



- 15.En utilisant l'outil de montage, placez la bague de rétention dans le logement de filtre et appuyez fermement pour la mettre en place.
- 16. Faites tourner l'outil jusqu'à ce que l'encoche de la bague de rétention soit alignée avec l'extrémité de l'outil de montage de filtre, puis retirez l'outil.



- 17.Replacez la roue porte-filtres dans son logement et enfoncez-la jusqu'à ce qu'elle soit fixée magnétiquement de manière sûre.
- 18.Replacez la plaque de protection sur le dessous de l'instrument en la fixant avec les trois vis orange.
- 19. Replacez l'appareil dans sa position normale.
- 20.Cliquez sur **OK** pour finaliser la procédure et initialiser la roue porte-filtres.
- 21.Définissez le filtre nouvellement inséré (veuillez consulter le chapitre suivant pour la procédure détaillée).



# 7.5.2 Définir les filtres

Dans la boîte de dialogue **Définition du filtre**, assignez les longueurs d'onde appropriées aux filtres remplacés en entrant les nouvelles longueurs d'onde aux endroits qui conviennent.

Si un filtre a été inséré à une nouvelle position, activez la position correspondante en sélectionnant la case à cocher et en entrant la longueur d'onde appropriée. En cliquant sur **Enregistrer**, les définitions de filtres seront enregistrées et les filtres seront initialisés.

Une fois les filtres initialisés, l'instrument est prêt pour les mesures.



Note Veillez à ne pas mélanger les positions des filtres et les longueurs d'onde de filtre, sous peine de fausser les données de mesure.

# 7.6 Élimination des déchets

# 7.6.1 Introduction

Suivez les procédures de laboratoire pour l'élimination des déchets présentant un danger biologique, conformément aux régulations locales et nationales.

Ce chapitre fournit des instructions expliquant comment éliminer le matériel de déchet produit en lien avec l'INFINITE F50 PLUS, dans le respect de la loi.



#### PRÉCAUTION

RESPECTEZ TOUTES LES RÉGULATIONS ENVIRONNEMENTALES NATIONALES, RÉGIONALES ET LOCALES.

# 7.6.2 Élimination du matériel d'emballage

Le matériel d'emballage est fait de matériaux recyclables. Si vous ne souhaitez pas le conserver en vue d'utilisations futures, par ex. à des fins de transport et de stockage, éliminez le matériel d'emballage conformément aux régulations locales.

# 7.6.3 Élimination du matériel de fonctionnement

#### AVERTISSEMENT

DES DANGERS BIOLOGIQUES PEUVENT DÉRIVER DU MATÉRIEL DE DÉCHET (MICROPLAQUES) DU PROCESSUS EXÉCUTÉ AVEC LE LECTEUR D'ABSORBANCE INFINITE F50 PLUS.



TRAITEZ LES MICROPLAQUES UTILISÉES, LES AUTRES CONSOMMABLES ET TOUTES LES SUBSTANCES EMPLOYÉES DANS LE RESPECT DES BONNES PRATIQUES DE LABORATOIRE.

RENSEIGNEZ-VOUS SUR LES LIEUX DE COLLECTE ET SUR LES MÉTHODES D'ÉLIMINATION APPROUVÉES DANS VOTRE PAYS OU VOTRE RÉGION.



## 7.6.4 Élimination de l'instrument

Si vous avez des questions concernant l'élimination de l'instrument, veuillez contacter votre représentant Tecan local.

Degré de pollution	2 (IEC/EN 61010-1)
Méthode d'élimination	Déchets contaminés





DIRECTIVE 2012/19/UE SUR LES DÉCHETS D'ÉQUIPEMENTS ÉLECTRIQUES ET ÉLECTRONIQUES (DEEE) EFFETS NÉGATIFS SUR L'ENVIRONNEMENT ASSOCIES AU TRAITEMENT DES DÉCHETS. • NE TRAITEZ PAS LES DÉCHETS D'ÉQUIPEMENTS ÉLECTRIQUES ET ÉLECTRONIQUES COMME DES DÉCHETS MÉNAGERS NON-TRIES.

ATTENTION

• COLLECTEZ SÉPARÉMENT LES DÉCHETS D'ÉQUIPEMENTS ÉLECTRIQUES ET ÉLECTRONIQUES.

#### AVERTISSEMENT

SELON LES APPLICATIONS, CERTAINES PARTIES DE L'INFINITE F50 PLUS PEUVENT AVOIR ÉTÉ EN CONTACT AVEC DES PRODUITS PRÉSENTANT UN RISQUE BIOLOGIQUE.

- ASSUREZ-VOUS DE TRAITER CE MATÉRIEL CONFORMÉMENT AUX STANDARDS ET AUX RÉGULATIONS DE SÉCURITÉ APPLICABLES.
- DÉCONTAMINEZ TOUTES LES PARTIES AVANT L'ÉLIMINATION.

Pour plus d'informations sur le produit, veuillez contacter Tecan Austria GmbH Untersbergstrasse 1A A-5082 Grödig/Salzbourg AUTRICHE/EUROPE T +43 6246 8933 444 E-mail : <u>expertline-at@tecan.com</u> www.tecan.com


# 8. Dépannage

## 8.1 Introduction

Le microprocesseur interne contrôle et vérifie toutes les fonctions électroniques, ainsi que les mesures, les opérations et les résultats. Si le microprocesseur détecte une erreur ou une procédure de fonctionnement incorrecte, un message d'erreur s'affiche sur l'ordinateur.

## 8.1.1 Tableau des messages d'erreur et de dépannage

Le tableau suivant fournit une brève description des messages d'erreur et des actions de dépannage.



Note Si des messages d'erreur non cités ci-dessous s'affichent, veuillez contacter votre représentant Tecan local.

Message d'erreur	Description	Dépannage
Erreur de système		
"Lid Open Error" (Erreur : couvercle ouvert)	Couvercle ouvert en début de mesure	Fermez le couvercle et redémarrez la mesure
"MTP Init Error" (Erreur d'initialisation de plaque MT)	Impossible d'initialiser le transport de plaque MT	Problème matériel : défaut électronique, courroie rompue ou blocage mécanique du transport de plaque MT
"MTP lost steps abs( <i>steploss</i> ) > max_steploss" (La plaque MT a perdu des pas abs(perte_pas) > perte_pas_max)	La plaque MT a perdu des pas au cours de la mesure <i>perte_pas</i> : nombre de pas perdus <i>perte_pas_max</i> : nombre de pas perdus autorisé	Problème matériel : défaut électronique, mécanique grippée
"Filter lost steps abs( <i>steploss</i> ) > <i>max_steploss</i> " (Le filtre a perdu des pas abs(perte_pas) > perte_pas_max)	La roue porte-filtres a perdu des pas pendant la mesure. <i>perte_pas</i> : nombre de pas perdus <i>perte_pas_max</i> : nombre de pas perdus autorisé	Problème matériel : Electronique déficiente, mécanique grippée
"USB timeout" (Dépassement de délai USB)	Dépassement de délai de la communication USB	Erreur système – consulter l'assistance client
"Lamp Low! Minimum: <i>minimum</i> , Maximum: <i>maximum</i> " (Lampe faible ! Minimum : minimum, maximum : maximum) <i>Ce message s'affiche avec les versions V1.11 et antérieures du</i> <i>micrologiciel.</i>	L'intensité de lumière mesurée n'atteint pas l'intervalle attendu situé entre <i>minimum</i> et <i>maximum</i>	Problème matériel : Défaut électronique, fibre cassée



### 8. Dépannage

Message d'erreur	Description	Dépannage
À partir de la version V1.12 du micrologiciel, le message "Lamp Low" est remplacé par : "Prepare REF check"	L'intensité de lumière mesurée n'atteint pas l'intervalle attendu situé entre <i>minimum</i> et <i>maximum</i>	Problème matériel : Défaut électronique, fibre cassée
"Wavelength Not Available ! Wavelength: <i>wavelength</i> nm" (Longueur d'onde indisponible ! Longueur d'onde : longueur d'onde nm)	Le filtre de la longueur d'onde <i>longueur d'onde</i> n'a pas été trouvé dans la roue porte-filtres	Erreur système – consulter l'assistance client
"Channel Low! Channel: <i>channel_nr</i> , Minimum: <i>minimum</i> , Maximum: <i>maximum</i> " (Canal faible ! Canal : canal_nr, minimum : minimum, maximum : maximum)	Le signal du canal <i>canal_nr</i> n'a pas atteint la plage située entre <i>minimum</i> et <i>maximum</i>	Problème matériel : défaut électronique, fibre cassée
"Invalid Wavelength! Wavelength: <i>wavelength</i> nm" (Longueur d'onde non valide ! Longueur d'onde : longueur d'onde nm)	La longueur d'onde du filtre est hors de l'intervalle de longueur d'onde des DEL blanche et bleue.	Erreur système – consulter l'assistance client
"Lamp Overflow!" Minimum: <i>minimum</i> , Maximum: <i>maximum</i> " (Débordement lampe ! Minimum : minimum, maximum : maximum)	Le signal sur l'ADC dépasse la plage attendue entre <i>minimum</i> et <i>maximum</i>	Problème matériel : défaut électronique
"Value Not Set: <i>value -1</i> " (Valeur non définie : valeur -1)	La valeur <i>valeur</i> n'est pas définie	Erreur système – consulter l'assistance client
"Filter Init Error" (Erreur d'initialisation de filtre)	Impossible d'initialiser le transport de filtre	Problème matériel : défaut électronique, transport de roue porte-filtres bloqué mécaniquement

## 8.1.2 Définition de 'Overflow' (débordement)

Si le résultat des mesures d'absorbance est en dehors des spécifications de l'instrument (> 4.0 DO), un débordement peut survenir, et la valeur de DO mesurée du puits en cours sera remplacée par 'Overflow'. Ceci est effectué par le logiciel de contrôle, et non par l'instrument lui-même.

## 8.1.3 Panne électrique

Voici ce qui arrive en cas de panne électrique :

Panne électrique de l'instrument, mais pas de l'ordinateur de contrôle (par ex. si l'ordinateur est branché à un onduleur) : la connexion USB entre l'instrument et l'ordinateur sera perdue. Un message d'erreur du logiciel de contrôle est créé. Panne électrique de l'instrument et de l'ordinateur de contrôle : l'ordinateur doit redémarrer. Aucune donnée de mesure ne sera disponible.



# 9. Abréviations, marques et symboles

## 9.1 Abréviations

Les abréviations suivantes sont fournies pour servir de référence, et peuvent apparaître dans la notice d'utilisation.

А	Ampère
AC	Courant alternatif
ADC	Convertisseur analogique digital
ANSI/SBS	American National Standards Institute/Society for Biomolecular Screening (Institut américain de normalisation/Société des sciences Biomoléculaires)
ASCII	American Standard Code for Information Interchange (Code américain normalisé pour l'échange d'information)
ASTM	American Society for Testing and Material (Société américaine pour les essais et les matériaux)
°C	Degrés Celsius
CE	Marquage CE de conformité
CFR	Code of Federal Regulations (Code des règlements fédéraux)
cm	Centimètre
DC	Courant continu
EC	Communauté européenne
ELISA	Dosage immunoenzymatique sur support solide
EN	Norme européenne
°F	Degrés Fahrenheit
FDA	Food and Drug Administration (Administration des médicaments des denrées alimentaires et des médicaments)
Hz	Hertz
IEC (CEI)	International Electrotechnical Commission (Commission électrotechnique internationale)
ID	Identification
IFU	Notice d'utilisation
IQ	Qualification de l'installation
IVD	Diagnostic in vitro
IVDR	Règlementation de diagnostic in vitro (RDIV) (UE) 2017/746
kg	Kilogramme
1	Litre
DEL	Diode électro-luminescente
LIS	Laboratory Information System (Système d'information de laboratoire)
mg	Milligramme
ml	Millilitre
mm	Millimètre



#### 9. Abréviations, marques et symboles

MTP	Microplaque
μΙ	Microlitre
MNF	Métal non ferreux
NIST	National Institute of Standards and Technology (Institut national des standards et de la technologie)
nm	nanomètre
NRTL	Nationally Recognized Testing Laboratory (Laboratoire de test reconnu au niveau national)
DO	Densité optique
QO	Qualification opérationnelle
PCB	Printed Circuit Board (circuit imprimé)
RF	Radio fréquence
RoHS	Restriction of the Use of Certain Hazardous Substances (Restriction de l'utilisation de certaines substances dangereuses)
POP	Procédures opérationnelles permanents
USB	Universal Serial Bus (Bus série universel)
UA	Unité arbitraire
TÜV	Technischer Überwachungsverein (Agence d'inspection technique)
V	Volt
VA	Volt ampère
WEEE	Waste electrical and electronic equipment (Déchets d'équipements électriques et électroniques)

## 9.2 Marques de commerce

Les noms de produits suivants et toutes les marques déposées et non déposées mentionnées dans ce document ne sont utilisés qu'à des fins d'identification et restent la propriété exclusive de leurs détenteurs respectifs.

- Magellan<sup>™</sup>, Infinite<sup>®</sup>, MultiCheck<sup>™</sup>, Tecan<sup>®</sup> et le logo Tecan sont des marques déposées de Tecan Group Ltd., Männedorf, Suisse
- Windows® et Excel® sont des marques déposées de Microsoft Corporation, Redmond, WA, USA
- Pentium<sup>®</sup> et Atom<sup>™</sup> sont des marques déposées d'Intel Corporation, Santa Clara, CA, USA
- Adobe® Reader® est une marque déposée d'Adobe Systems Incorporated, Seattle, WA, USA
- Microcide SQ<sup>™</sup> est une marque commerciale de Global Biotechnologies Inc., Portland, ME, USA
- Decon 90<sup>™</sup> est une marque commerciale de Decon Laboratories Ltd., Hove, East Sussex, Grande-Bretagne



## 9.3 Symboles

Les symboles suivants apparaissent dans ce document.

	Fabricant
M	Date de fabrication
CE	Marquage CE de conformité
UK CA	United Kingdom Conformity Assessed Le marquage United Kingdom Conformity Assessed indique la conformité du produit étiqueté aux exigences du règlement applicable en Grande-Bretagne.
Ĩ	Consulter le manuel de l'utilisateur
IVD	Dispositif médical de diagnostic in vitro
UDI	Unique Device Identification – Identifiant unique des dispositifs Le symbole UDI identifie le support de données sur l'étiquette.
REF	Numéro da catalogue
SN	Numéro de série
€ ↓	Symbole USB
X	Symbole WEEE
	Symbole RoHS, Chine
	NRTL TÜV SÜD MARK
	Risques biologiques



# Index

## Α

Abréviations	75
Absorbance	23
Agitation	23, 33
Assistant	
Attacher une signature	19
Créer/éditer une liste ID d'échantillon	19
Créer/éditer une méthode	19
Démarrer la mesure	19
Liste des assistants	19
Assistant Evaluer les résultats	19
Attendre (horloge)	25

### В

2
23
24
25
23
22
23
24
25
23
22

#### С

Certificat de sécurité	66
Changement de filtre	68
Commentaire	24
Condition cinétique	24
Contrôle qualité	39
Cycle cinétique	24

## D

Déballage	
déballage et inspection	
procédure de déballage	
Débordement	74
Définir les filtres	71
Définir les mesures	
Dépannage	73
Désinfection	63, 64
Certificat de sécurité	
Procédure	65
Solutions	64

## Ε

Editeur de paramètres de mesure	21
Elements de programme	
Indentation et libération	30
Partie de la plaque	22
Eléments de programme	
Absorbance	23
Agitation	23

Attendre (horloge)	25
Commentaire	24
Condition cinétique	24
Cycle cinétique	24
Incubation	25
Plaque	22
Requête de l'utilisateur	25
Elimination	
Matériel de fonctionnement	71
Élimination	
Instrument	72
Matériel d'emballage	71
Elimination des déchets	63
Élimination des déchets	
Exemple d'application	43
ELISA quantitatif	
- 1	

## I

Incubation	
Instrument	
Accessoires	
Description	33
Emplacement	32
Fonctionnalités	33
Spécifications	

## L

Liste des assistants	19
Logiciel	17
Installation	17
Interface utilisateur	19
Liste des assistants	19
Magellan	17
Magellan Tracker	17
Qualification de l'installation	18

## М

17
17
63, 67
67
73
29
27
28
16
38
32

### Ν

Nécessités d'alimentation	14
Nécessités environnementales	14
Nettoyage	63

### 0

Optimisation des performances ...... 32



## Index

Ρ

Panne électrique	
Partie de la plaque	
Plaque	
Procédure d'auto-vérification	
Profils d'utilisateur	11
-	

### Q

Qualification opérationnelle	39
Test de linéarité	41
Test de microplaque	39
Test de précision	40
Test MultiCheck	39

## R

Remplacement du filtre Requête de l'utilisateur Roue porte-filtres	67 25 35
S	
Sécurité Spécifications Symboles Système nécessaire	7 36 77 15
V	
Volet d'info	26



# **Service client Tecan**

Si vous avez des questions ou si vous avez besoin d'une assistance technique pour votre produit Tecan, veuillez contacter votre service client Tecan local. Rendez-vous à l'adresse <u>http://www.tecan.com/</u> pour obtenir les coordonnées de contact.

Avant de contacter Tecan, préparez les informations suivantes afin d'obtenir le meilleur support technique possible (voir la plaque signalétique) :

- Nom de modèle de votre produit
- Numéro de série (SN) de votre produit
- Logiciel et version logicielle (le cas échéant)
- Description du problème et interlocuteur
- Date et heure de la survenue du problème
- Mesures déjà prises pour corriger le problème
- Vos coordonnées de contact (numéro de téléphone, numéro de fax, adresse 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.