Οδηγίες χρήσης για το **INFINITE F50 PLUS**

μετρητή απορρόφησης

MAGELLAN

λογισμικό

•TECAN•









ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΔΙΑΒΑΣΤΕ ΠΡΟΣΕΚΤΙΚΑ ΚΑΙ ΑΚΟΛΟΥΘΗΣΤΕ ΤΙΣ ΟΔΗΓΙΕΣ ΧΡΗΣΗΣ ΠΡΙΝ ΤΗ ΛΕΙΤΟΥΡΓΙΑ ΤΟΥ ΟΡΓΑΝΟΥ.

Σημείωση

Έχει καταβληθεί κάθε δυνατή προσπάθεια για την αποφυγή σφαλμάτων στο κείμενο και τα διαγράμματα. Ωστόσο, η Tecan Austria GmbH δεν αναλαμβάνει καμία ευθύνη για τυχόν σφάλματα στο παρόν έγγραφο.

Πολιτική της Tecan Austria GmbH αποτελεί η βελτίωση των προϊόντων, καθώς νέες τεχνικές και στοιχεία γίνονται διαθέσιμα. Κατά συνέπεια, η Tecan Austria GmbH διατηρεί το δικαίωμα της αλλαγής των προδιαγραφών οποιαδήποτε στιγμή με την κατάλληλη επαλήθευση, επικύρωση και εγκρίσεις.

Εκτιμούμε την παροχή σχολίων εκ μέρους σας σχετικά με το παρόν έγγραφο.



Κατασκευαστής

Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria T: +43 6246 89330

www.tecan.com E-mail: office.austria@tecan.com

Πληροφορίες πνευματικών δικαιωμάτων

Τα περιεχόμενα του παρόντος εγγράφου αποτελούν ιδιοκτησία της Tecan Austria GmbH. Απαγορεύεται η αντιγραφή, η αναπαραγωγή ή η μεταβίβαση σε άλλα πρόσωπα χωρίς προηγούμενη έγγραφη άδεια.

Copyright © Tecan Austria GmbH Με την επιφύλαξη παντός δικαιώματος. Εκτυπώθηκε στην Αυστρία.

Δήλωση για Πιστοποίηση Ε.Ε.

Ανατρέξτε στην τελευταία σελίδα του παρόντος.

Πληροφορίες σχετικά με τις Οδηγίες Χρήσης.

Πρωτότυπες οδηγίες. Το παρόν έγγραφο προορίζεται να χρησιμοποιηθεί ως **Οδηγίες Χρήσης** (IFU) για το μετρητή απορρόφησης INFINITE F50 PLUS, ο οποίος έχει σχεδιαστεί για τη μέτρηση της απορρόφησης φωτός (οπτικής πυκνότητας) δειγμάτων σε μικροπλάκες 96 κυψελίδων. Προορίζεται να χρησιμοποιηθεί ως αναφορά και οδηγίες για το χρήστη.

Στο έγγραφο παρέχονται οδηγίες για:

- Εγκατάσταση του οργάνου
- Λειτουργία του οργάνου
- Καθαρισμό και συντήρηση του οργάνου

Παρατηρήσεις για τις εικόνες από το λογισμικό

Ο αριθμός έκδοσης που αναφέρεται στις εικόνες ενδέχεται να μην ανταποκρίνεται πάντα στην πιο πρόσφατη έκδοση. Οι εικόνες αντικαθίστανται μόνο εάν το περιεχόμενο που σχετίζεται με την εφαρμογή έχει αλλάξει.



Ιστορικό αλλαγών – Οδηγίες χρήσης

Έκδοση	Αλλαγές
V 1.2	Νέα Δήλωση συμμόρφωσης (DoC)

Επισημάνσεις προειδοποίησης, προσοχής και σημειώσεων

Οι παρακάτω επισημάνσεις χρησιμοποιούνται σε όλο το παρόν έγγραφο για να υποδείξουν σημαντικές πληροφορίες ή να προειδοποιήσουν το χρήστη για πιθανώς επικίνδυνες καταστάσεις:

> Σημείωση Παρέχει χρήσιμες πληροφορίες.

ΠΡΟΣΟΧΗ ΥΠΟΔΕΙΚΝΥΕΙ ΤΗΝ ΠΙΘΑΝΟΤΗΤΑ ΒΛΑΒΗΣ ΣΤΟ ΟΡΓΑΝΟ Ή ΑΠΩΛΕΙΑΣ ΔΕΔΟΜΕΝΩΝ ΣΕ ΠΕΡΙΠΤΩΣΗ ΠΟΥ ΔΕΝ ΤΗΡΗΘΟΥΝ ΟΙ ΟΔΗΓΙΕΣ.

ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΥΠΟΔΕΙΚΝΥΕΙ ΤΗΝ ΠΙΘΑΝΟΤΗΤΑ ΣΟΒΑΡΟΥ ΠΡΟΣΩΠΙΚΟΥ ΤΡΑΥΜΑΤΙΣΜΟΥ, ΑΠΩΛΕΙΑΣ ΖΩΗΣ Ή ΒΛΑΒΗΣ ΣΤΟΝ ΕΞΟΠΛΙΣΜΟ ΣΕ ΠΕΡΙΠΤΩΣΗ ΠΟΥ ΔΕΝ ΤΗΡΗΘΟΥΝ ΟΙ ΟΔΗΓΙΕΣ.

ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΥΠΟΔΕΙΚΝΥΕΙ ΤΗΝ ΠΙΘΑΝΗ ΠΑΡΟΥΣΙΑ ΒΙΟΛΟΓΙΚΑ ΕΠΙΚΙΝΔΥΝΟΥ

ΑΣΦΑΛΕΙΑΣ ΕΡΓΑΣΤΗΡΙΟΥ.















ΠΡΟΣΟΧΗ

ΟΔΗΓΙΑ 2012/19/ΕΕ ΠΕΡΙ ΑΠΟΒΛΗΤΩΝ ΗΛΕΚΤΡΙΚΟΥ ΚΑΙ ΗΛΕΚΤΡΟΝΙΚΟΥ ΕΞΟΠΛΙΣΜΟΥ (WEEE)

ΑΡΝΗΤΙΚΕΣ ΠΕΡΙΒΑΛΛΟΝΤΙΚΕΣ ΕΠΙΠΤΩΣΕΙΣ ΠΟΥ ΣΧΕΤΙΖΟΝΤΑΙ ΜΕ ΤΗ ΔΙΑΧΕΙΡΙΣΗ ΑΠΟΒΛΗΤΩΝ.

- ΜΗΝ ΔΙΑΧΕΙΡΙΖΕΣΤΕ ΗΛΕΚΤΡΙΚΟ ΚΑΙ ΗΛΕΚΤΡΟΝΙΚΟ ΕΞΟΠΛΙΣΜΟ ΩΣ ΟΙΚΙΑΚΑ ΑΠΟΒΛΗΤΑ.
- ΣΥΛΛΕΓΕΤΕ ΞΕΧΩΡΙΣΤΑ ΤΑ ΑΠΟΒΛΗΤΑ ΑΠΟ ΗΛΕΚΤΡΙΚΟ ΚΑΙ ΗΛΕΚΤΡΟΝΙΚΟ ΕΞΟΠΛΙΣΜΟ.



Πίνακας περιεχομένων

1.	Ασφ	αλεια		7
	1.1	Ασφάλ	ιεια οργάνου	7
2.	Γενι	κά		g
_ .	2 1	Поосо	νζόμενη νοήση	
	2.1	Λονή Λ	ηςομενή χρηση	
	2.2	πρχη /	λιορστών	10
	2.5	224	Έμπειοος γρήστης Επίπεδο διαγειοιστή	
		2.3.1	Εμπειρος χρηστης – Εππεσο σιαχειριστη	
		2.3.2	Τελικός χρηστης η χρηστης «ροστίνας»	
•	-	2.3.3	τεχνικός σερρις	
3.	Ιατ	τρωτα β	ηματα	
	3.1	Άνοιγμ	ια συσκευασίας και επιθεώρηση	
		3.1.1	Επιθεώρηση της συσκευασίας που σας παραδόθηκε	
		3.1.2	Διαδικασία ανοίγματος συσκευασίας	13
	3.2	Απαιτή	ήσεις τροφοδοσίας	14
	3.3	Περιβα	ιλλοντικές απαιτήσεις	14
	3.4	Απαιτή	ήσεις συστήματος	15
	3.5	Ενεργα	οποίηση του οργάνου	16
	3.6	Λογισμ	JIKÓ	17
		3.6.1	Εισαγωγή / Προοριζόμενη χρήση	17
		3.6.2	Διαδικασία εγκατάστασης	17
		3.6.3	Επαλήθευση εγκατάστασης	
		3.6.4	Τα πρώτα βήματα με το Magellan	
	3.7	Magell	an – Πρόγραμμα επεξεργασίας παραμέτρων μέτρησης	21
		3.7.1	Γραμμή ελέγχου	
		3.7.2	Παράθυρο ροής εργασιών	
		3.7.3	Παράθυρο πληροφοριών	27
	3.8	Magell	an – Ορισμός μετρήσεων	
		3.8.1	Ορισμός μετρήσεων τελικού σημείου	
		3.8.2	Ορισμός μετρήσεων πολλαπλής σήμανσης	
		3.8.3	Ορισμός μετρήσεων ανάλυσης κινητικής	
		3.8.4	Δέσμευση και αποδέσμευση στοιχείων προγράμματος	
	3.9	Βελτισ	τοποίηση για καλύτερη απόδοση	
		3.9.1	Θέση οργάνου	
		3.9.2	Διαδικασία λειτουργίας	
4.	Λειτ	ουργίες	οργάνου	35
	4.1	Λειτου	ργίες οργάνου	35
		4.1.1	Ανακίνηση μικροπλάκας	35
	4.2	Περιγρ	αφή οργάνου	35
	4.3	Περιγρ	αφή τροχού φίλτρων	37
	4.4	Προδια	αγραφές οργάνου	
		4.4.1	Γενικές προδιαγραφές	
		4.4.2	Προδιαγραφές μέτρησης	
		4.4.3	Μικροπλάκες	
	4.5	Παρελι	κόμενα οργάνου	40
5.	Έλε	γχος πο	ριότητας	41
	5.1	Εισαγά	υγή	41
	5.2	Διαδικ	ασία αυτόματου ελέγχου	41
	5.3	Επαλή	θευση λειτουργικότητας	41
		5.3.1	Έλεγχος πολλαπλών σημείων (MultiCheck)	41
		5.3.2	Έλεγχος μικροπλάκας	41
		5.3.3	Ακρίβεια οργάνου με υγρά δείγματα	42
		5.3.4	Γραμμικότητα οργάνου με υγρά δείγματα	43
6.	Παο	άδεινυα	εφαρμογής	45
	6.1	Εισανι	۵۷ή	
	6.2	Παράδ	εινμα «βήμα-προς-βήμα»: Ποσοτικός προσδιορισμός ELISA	45



		6.2.1	Περιγραφή κιτ ελέγχου	
		6.2.2	Δημιουργία μεθόδου	
		6.2.3	Οργάνωση έντυπης αναφοράς	
		6.2.4	Εκτέλεση της μεθόδου	
		6.2.5	Αξιολόγηση αποτελέσματος	
		6.2.6	Σύνοψη ορισμού ποσοτικού προσδιορισμού ELISA στο Magellan	
7.	Καθα	αρισμός	, συντήρηση και απόρριψη	67
	7.1	Εισαγω	υγή	
	7.2	Καθαρι	ισμός του οργάνου	
	7.3	Απολύ	μανση οργάνου	
		7.3.1	Διαλύματα απολύμανσης	
		7.3.2	Διαδικασία απολύμανσης	
		7.3.3	Πιστοποιητικό ασφάλειας	
	7.4	Πρόγρα	αμμα προληπτικής συντήρησης για το INFINITE F50 PLUS	71
		7.4.1	Μηνιαία	71
		7.4.2	Κάθε 4 έτη	71
	7.5	Αντικατ	τάσταση και τοποθέτηση φίλτρων	71
		7.5.1	Διαδικασία αλλαγής φίλτρων	
		7.5.2	Ορισμός φίλτρων	75
	7.6	Απόρρ	ıψη	
		7.6.1	Εισαγωγή	
		7.6.2	Απόρριψη υλικών συσκευασίας	
		7.6.3	Απόρριψη υλικών λειτουργίας	
		7.6.4	Απόρριψη του οργάνου	
8.	Επίλ	υση πρ	οβλημάτων	77
	8.1	Εισαγω	υνή	
		8.1.1	Πίνακας μηνυμάτων σφάλματος και επίλυση προβλημάτων	
		8.1.2	Ορισμός «υπερχείλισης»	
		8.1.3	Διακοπή τροφοδοσίας	
9.	Συντ	ομογρα	φίες, εμπορικά σήματα και σύμβολα	79
	9.1	 Συντομ	ονραφίες	
	9.2	Εμπορ	ικά σήματα	
	9.3	Σύμβο λ	 \α	
Ευρε	τήριο	· ·		83
FEum	ποέτη	ιση πελ	ατών της Tecan	85
->	IPCII			



1. Ασφάλεια

1.1 Ασφάλεια οργάνου

- Τηρείτε πάντα τα βασικά μέτρα ασφαλείας κατά τη χρήση του προϊόντος ώστε να μειώσετε τον κίνδυνο τραυματισμού, πυρκαγιάς ή ηλεκτροπληξίας.
- Διαβάστε και βεβαιωθείτε ότι έχετε κατανοήσει όλες τις πληροφορίες στις Οδηγίες Χρήσης (IFU).
 Η μη ανάγνωση, κατανόηση και τήρηση των οδηγιών του παρόντος εγγράφου

Η μη αναγνωση, κατανοηση και τηρηση των σοηγιων του παροντος εγγραφου ενδέχεται να έχουν ως αποτέλεσμα βλάβη στο προϊόν, τραυματισμό του προσωπικού που το χειρίζεται ή χαμηλή απόδοση του οργάνου. Η Tecan δεν είναι υπεύθυνη για βλάβη ή τραυματισμούς που προέρχονται από λανθασμένο χειρισμό της συσκευής.

- Τηρείτε όλες τις επισημάνσεις με την ένδειξη ΠΡΟΕΙΔΟΠΟΙΗΣΗ και ΠΡΟΣΟΧΗ του παρόντος εγγράφου.
- Αποσυνδέετε πάντα τη συσκευή από την κύρια παροχή τροφοδοσίας πριν τον καθαρισμό και την απολύμανση.
- 5. Μην ανοίγετε το περίβλημα του οργάνου.
- 6. Τηρείτε όλα τα απαραίτητα μέτρα ασφαλείας εργαστηρίου, όπως η χρήση προστατευτικού ρουχισμού (π.χ. γάντια, μπλούζα εργαστηρίου και προστατευτικά γυαλιά) και εφαρμόζετε όλες τις εγκεκριμένες διαδικασίες ασφαλείας εργαστηρίου.



ΠΡΟΣΟΧΗ

ΕΑΝ ΟΙ ΟΔΗΓΙΕΣ ΠΟΥ ΠΕΡΙΕΧΟΝΤΑΙ ΣΤΟ ΠΑΡΟΝ ΕΓΓΡΑΦΟ ΔΕΝ ΤΗΡΗΘΟΥΝ, ΕΝΔΕΧΕΤΑΙ ΝΑ ΠΡΟΚΛΗΘΕΙ ΒΛΑΒΗ ΣΤΟ ΟΡΓΑΝΟ Ή ΟΙ ΔΙΑΔΙΚΑΣΙΕΣ ΝΑ ΜΗΝ ΕΚΤΕΛΕΣΤΟΥΝ ΣΩΣΤΑ ΚΑΙ ΔΕΝ ΜΠΟΡΟΥΜΕ ΝΑ ΕΓΓΥΗΘΟΥΜΕ ΤΗΝ ΑΣΦΑΛΕΙΑ ΤΟΥ ΟΡΓΑΝΟΥ.

Θεωρείται δεδομένο ότι οι χειριστές του οργάνου, εξαιτίας της επαγγελματικής τους εμπειρίας, είναι εξοικειωμένοι με τα απαραίτητα μέτρα ασφαλείας για το χειρισμό χημικών και βιολογικά επικίνδυνων ουσιών.

Τηρείτε την παρακάτω νομοθεσία και οδηγίες:

- Εθνική νομοθεσία για βιομηχανική προστασία
- Κανονισμούς αποφυγής ατυχημάτων
- Δελτία δεδομένων ασφαλείας των κατασκευαστών αντιδραστηρίων







ΑΝΑΛΟΓΑ ΜΕ ΤΙΣ ΕΦΑΡΜΟΓΕΣ, ΕΞΑΡΤΗΜΑΤΑ ΤΟΥ INFINITE F50 PLUS ΕΝΔΕΧΕΤΑΙ ΝΑ ΕΡΘΟΥΝ ΣΕ ΕΠΑΦΗ ΜΕ ΒΙΟΛΟΓΙΚΑ ΕΠΙΚΙΝΔΥΝΟ/ΜΟΛΥΣΜΑΤΙΚΟ ΥΛΙΚΟ.

ΒΕΒΑΙΩΘΕΙΤΕ ΟΤΙ ΤΟ ΟΡΓΑΝΟ ΧΡΗΣΙΜΟΠΟΙΕΙΤΑΙ ΜΟΝΟ ΑΠΟ ΚΑΤΑΛΛΗΛΑ ΕΚΠΑΙΔΕΥΜΕΝΟ ΠΡΟΣΩΠΙΚΟ. ΣΕ ΠΕΡΙΠΤΩΣΗ ΣΥΝΤΗΡΗΣΗΣ/ΕΠΙΣΚΕΥΗΣ Ή ΑΛΛΑΓΗΣ ΘΕΣΗΣ Ή ΑΠΟΡΡΙΨΗΣ ΤΟΥ ΟΡΓΑΝΟΥ, ΑΠΟΛΥΜΑΙΝΕΤΕ ΠΑΝΤΑ ΤΟ ΟΡΓΑΝΟ ΣΥΜΦΩΝΑ ΜΕ ΤΙΣ ΟΔΗΓΙΕΣ ΠΟΥ ΔΙΝΟΝΤΑΙ ΣΤΟ ΠΑΡΟΝ ΕΓΓΡΑΦΟ.

ΤΗΡΕΙΤΕ ΤΑ ΚΑΤΑΛΛΗΛΑ ΜΕΤΡΑ ΑΣΦΑΛΕΙΑΣ ΕΡΓΑΣΤΗΡΙΟΥ, ΟΠΩΣ Η ΧΡΗΣΗ ΠΡΟΣΤΑΤΕΥΤΙΚΟΥ ΡΟΥΧΙΣΜΟΥ ΚΑΤΑ ΤΗΝ ΕΡΓΑΣΙΑ ΜΕ ΠΙΘΑΝΩΣ ΜΟΛΥΣΜΑΤΙΚΕΣ ΟΥΣΙΕΣ.

ΠΡΟΕΙΔΟΠΟΙΗΣΗ

ΤΟ ΟΡΓΑΝΟ ΣΥΜΜΟΡΦΩΝΕΤΑΙ ΜΕ ΤΙΣ ΑΠΑΙΤΗΣΕΙΣ ΕΚΠΟΜΠΩΝ ΚΑΙ ΑΤΡΩΣΙΑΣ ΠΟΥ ΠΕΡΙΓΡΑΦΟΝΤΑΙ ΣΤΟ ΠΡΟΤΥΠΟ ΙΕС 61326-2-6. ΩΣΤΟΣΟ, ΤΟ ΗΛΕΚΤΡΟΜΑΓΝΗΤΙΚΟ ΠΕΡΙΒΑΛΛΟΝ ΘΑ ΠΡΕΠΕΙ ΝΑ ΑΞΙΟΛΟΓΗΘΕΙ ΠΡΙΝ ΤΗ ΛΕΙΤΟΥΡΓΙΑ ΤΟΥ ΟΡΓΑΝΟΥ.

Η ΕΞΑΣΦΑΛΙΣΗ ΤΗΡΗΣΗΣ ΣΥΜΒΑΤΟΥ ΗΛΕΚΤΡΟΜΑΓΝΗΤΙΚΟΥ ΠΕΡΙΒΑΛΛΟΝΤΟΣ ΩΣΤΕ ΤΟ ΟΡΓΑΝΟ ΝΑ ΛΕΙΤΟΥΡΓΕΙ ΣΥΜΦΩΝΑ ΜΕ ΤΗΝ ΠΡΟΟΡΙΖΟΜΕΝΗ ΧΡΗΣΗ ΑΠΟΤΕΛΕΙ ΕΥΘΥΝΗ ΤΟΥ ΧΕΙΡΙΣΤΗ.

ΜΗ ΧΡΗΣΙΜΟΠΟΙΕΙΤΕ ΤΟ ΟΡΓΑΝΟ ΚΟΝΤΑ ΣΕ ΠΗΓΕΣ ΙΣΧΥΡΗΣ ΗΛΕΚΤΡΟΜΑΓΝΗΤΙΚΗΣ ΑΚΤΙΝΟΒΟΛΙΑΣ (Π.Χ. ΣΚΟΠΙΜΕΣ ΠΗΓΕΣ ΡΑΔΙΟΣΥΧΝΟΤΗΤΑΣ ΧΩΡΙΣ ΘΩΡΑΚΙΣΗ) ΚΑΘΩΣ ΑΥΤΟ ΕΝΔΕΧΕΤΑΙ ΝΑ ΕΠΗΡΕΑΣΕΙ ΤΗ ΣΩΣΤΗ ΛΕΙΤΟΥΡΓΙΑ ΤΟΥ ΟΡΓΑΝΟΥ ΚΑΙ ΝΑ ΟΔΗΓΗΣΕΙ ΣΕ ΛΑΝΘΑΣΜΕΝΑ ΑΠΟΤΕΛΕΣΜΑΤΑ.



2. Γενικά

2.1 Προοριζόμενη χρήση

To INFINITE F50 PLUS είναι ένας αυτοματοποιημένος μετρητής απορρόφησης 96 κυψελίδων που περιλαμβάνει το λογισμικό Magellan για επαγγελματική χρήση σε εργαστήριο για τη μέτρηση της απορρόφησης φωτός (οπτική πυκνότητα) ομοιογενών υγρών μέσων σε in vitro διαγνωστική χρήση.

Το όργανο προορίζεται για χρήση κυρίως σε in vitro διαγνωστική ανάλυση δειγμάτων του ανθρώπινου σώματος με επιλεγμένες ενζυμικές ανοσοπροσροφητικες μεθόδους προσδιορισμού (ELISA). Οι συγκεκριμένες διαγνωστικές πληροφορίες και ο τύπος του δείγματος καθορίζονται από την επιλεγμένη ανάλυση.

The INFINITE F50 PLUS προορίζεται για τη μέτρηση και για την αξιολόγηση ποιοτικών και ποσοτικών, ημι-ποσοτικών, και ποσοτικών αναλύσεων σύμφωνα με τις προγραμματισμένες παραμέτρους διαγνωστικού ελέγχου και τις προδιαγραφές του οργάνου.

Το προϊόν προορίζεται για επαγγελματική εργαστηριακή χρήση από εξιδεικευμένο προσωπικό. Το προϊόν δεν προορίζεται για οικιακή χρήση ή χρήση από απλούς εργαζόμενους.

Σημείωση Εάν ο μετρητής απορρόφησης INFINITE F50 PLUS ή το λογισμικό Magellan τροποποιηθούν με οποιονδήποτε τρόπο, η εγγύηση δεν θα ισχύει και το όργανο δεν θα συμμορφώνεται με τους κανονισμούς.







Εάν ο μετρητής απορρόφησης INFINITE F50 PLUS χρησιμοποιηθεί διαφορετικά από ό,τι αναφέρεται στην ενότητα «Προοριζόμενη χρήση» παραπάνω ή χρησιμοποιηθεί με άλλο λογισμικό εκτός του Magellan, το όργανο δεν θα συμμορφώνεται πλέον με την οδηγία IVD και ο χρήστης θα είναι υπεύθυνος για την αντίστοιχη χρήση και την απαραίτητη επικύρωση.



Τα αποτελέσματα που λαμβάνονται με τη χρήση του INFINITE F50 PLUS επηρεάζονται από τη σωστή χρήση του οργάνου και των μικροπλακών, σύμφωνα με τις οδηγίες που παρέχονται στο παρόν έγγραφο, καθώς και από τις υγρές ενώσεις που χρησιμοποιούνται (αντιδραστήρια, χημικές αντιδράσεις). Οι οδηγίες χρήσης, αποθήκευσης και εφαρμογών που αφορούν δείγματα ή αντιδραστήρια πρέπει να ακολουθούνται αυστηρά. Κατά συνέπεια, τα αποτελέσματα θα πρέπει να ερμηνεύονται προσεκτικά.





Σημείωση Μην ανοίγετε το περίβλημα του οργάνου. Σε αυτήν την περίπτωση, η εγγύηση θα καταστεί άκυρη και δεν θα ισχύει.

2.2 Αρχή Λειτουργίας

To INFINITE F50 PLUS είναι ένας μετρητής απορρόφησης 8 καναλιών για τη μέτρηση της οπτικής πυκνότητας (OD) υγρών σε μικροπλάκες 96 κυψελίδων.

Το φως που δημιουργείται από τη λυχνία LED περνά μέσα από ένα οπτικό φίλτρο στον τροχό του φίλτρου για την επιλογή του μήκους κύματος που απαιτείται για τη μέτρηση της απορρόφησης.

Αφού περάσει μέσα από το φίλτρο, η δέσμη φωτός χωρίζεται σε οκτώ οπτικές ίνες για εστίαση σε οκτώ κυψελίδες της μικροπλάκας. Οι δέσμες φωτός που περνούν από τις κυψελίδες της μικροπλάκας, εστιάζονται από οκτώ οπτικούς φακούς και στη συνέχεια αποστέλλονται σε οκτώ φωτοδίοδους για ανίχνευση.

Αυτό σημαίνει ότι ο μετρητής INFINITE F50 PLUS μετρά οκτώ κυψελίδες παράλληλα, που αντιστοιχούν σε μία στήλη μιας μικροπλάκας 96 κυψελίδων.

Για τον έλεγχο και τη ρύθμιση της έντασης φωτός της λυχνίας LED, χρησιμοποιείται μία επιπλέον οπτική ίνα (κανάλι αναφοράς), όπου το φως παρακάμπτει τη μικροπλάκα.

Οι ληφθείσες τιμές διαπερατότητας μετατρέπονται σε τιμές οπτικής πυκνότητας (OD), σύμφωνα με τον ακόλουθο τύπο:

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

- I₀ = ένταση του προσπίπτοντος φωτός
- I = ένταση του διερχόμενου φωτός

Η τιμή OD είναι ο λογάριθμος του αντίστροφου της διαπερατότητας.

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



2.3 Προφίλ χρηστών

2.3.1 Έμπειρος χρήστης – Επίπεδο διαχειριστή

Ο διαχειριστής είναι αυτός που έχει την κατάλληλη τεχνική εκπαίδευση και τις αντίστοιχες δεξιότητες και εμπειρία. Εάν το προϊόν χρησιμοποιείται για τη χρήση για την οποία προορίζεται, το πρόσωπο αυτό μπορεί να αναγνωρίζει και να αποφεύγει κινδύνους.

Ο διαχειριστής διαθέτει εκτεταμένες δεξιότητες και μπορεί να καθοδηγήσει τον τελικό χρήστη ή το χρήστη «ρουτίνας» σε πρωτόκολλα προσδιορισμού σε σχέση με ένα προϊόν της Tecan και μέσα στα όρια τις προοριζόμενης χρήσης.

Απαιτούνται δεξιότητες χρήσης υπολογιστών και αγγλικής γλώσσας σε καλό επίπεδο.

2.3.2 Τελικός χρήστης ή χρήστης «ρουτίνας»

Ο τελικός χρήστης ή χρήστης «ρουτίνας» είναι αυτός που έχει την κατάλληλη τεχνική εκπαίδευση και τις αντίστοιχες δεξιότητες και εμπειρία. Εάν το προϊόν χρησιμοποιείται για τη χρήση για την οποία προορίζεται, το πρόσωπο αυτό μπορεί να αναγνωρίζει και να αποφεύγει κινδύνους.

Απαιτούνται δεξιότητες χρήσης εφαρμογών υπολογιστή και χειρισμού της γλώσσας για την αντίστοιχη τοπική γλώσσα στο χώρο εγκατάστασης, καθώς και γνώσεις Αγγλικών.

2.3.3 Τεχνικός σέρβις

Ο τεχνικός σέρβις είναι ένα πρόσωπο με την κατάλληλη τεχνική εκπαίδευση και τις αντίστοιχες δεξιότητες και εμπειρία. Εάν το προϊόν πρέπει να υποβληθεί σε σέρβις ή εργασίες συντήρησης, το πρόσωπο αυτό μπορεί να αναγνωρίζει και να αποφεύγει κινδύνους.

Απαιτούνται δεξιότητες χρήσης υπολογιστών και αγγλικής γλώσσας σε καλό επίπεδο.



Σημείωση Το τμήμα εξυπηρέτησης πελατών μπορεί να σας ενημερώσει για τις ημερομηνίες σεμιναρίων, καθώς και για τη διάρκεια και τη συχνότητά τους.

Διεύθυνση και αριθμό τηλεφώνου μπορείτε να βρείτε στο Internet: http://www.tecan.com/customersupport



3. Τα πρώτα βήματα

3.1 Άνοιγμα συσκευασίας και επιθεώρηση

3.1.1 Επιθεώρηση της συσκευασίας που σας παραδόθηκε

Το όργανο που σας παραδόθηκε περιλαμβάνει:

- Εξωτερικό τροφοδοτικό
- Καλώδιο τροφοδοσίας
- Καλώδιο USB για σύνδεση σε εξωτερικό υπολογιστή
- Οδηγίες Χρήσης, αρχεία PDF σε USB stick
- USB stick
 - Λογισμικό Magellan
 - Οδηγίες Χρήσης σε αρχεία PDF
 - ο Εργαλεία (π.χ. Adobe Reader)





Για να αποφύγετε ανεπιθύμητη απώλεια δεδομένων ή επίθεση από ιούς/κακόβουλο λογισμικό, μην καταργείτε την προστασία εγγραφής από το USB stick.

Σημείωση

ΠΡΟΣΟΧΗ

Η ΛΕΙΤΟΥΡΓΙΑ ΤΟΥ ΜΕΤΡΗΤΗ ΕΧΕΙ ΕΛΕΓΧΘΕΙ ΜΕ ΤΟ ΠΑΡΕΧΟΜΕΝΟ ΚΑΛΩΔΙΟ USB ΕΑΝ ΧΡΗΣΙΜΟΠΟΙΗΘΕΙ ΑΛΛΟ ΚΑΛΩΔΙΟ USB, ΔΕΝ ΜΠΟΡΟΥΜΕ ΝΑ ΕΓΓΥΗΘΟΥΜΕ ΤΗ ΣΩΣΤΗ ΑΠΟΔΟΣΗ ΤΟΥ ΟΡΓΑΝΟΥ

3.1.2 Διαδικασία ανοίγματος συσκευασίας

- 1. Πριν ανοίξετε τη συσκευασία, επιθεωρήστε την οπτικά για τυχόν ζημιά. Αναφέρετε αμέσως οποιαδήποτε ζημιά.
- 2. Επιλέξτε μια θέση για να τοποθετήσετε το όργανο. Η θέση θα πρέπει να είναι επίπεδη, χωρίς κραδασμούς, μακριά από άμεση επαφή με το ηλιακό φως και χωρίς σκόνη, διαλύτες και όξινους ατμούς. Βεβαιωθείτε ότι η απόσταση μεταξύ του οργάνου και του τοίχου ή άλλου εξοπλισμού είναι τουλάχιστον 5 cm.
- Ανασηκώστε το όργανο μέσα από τη συσκευασία και τοποθετήστε το στην επιλεγμένη θέση. Ανασηκώστε το όργανο με προσοχή.
- Επιθεωρήστε οπτικά το όργανο για εξαρτήματα που έχουν χαλαρώσει, λυγίσει ή σπάσει.
 Αναφέρετε αμέσως οποιαδήποτε ζημιά.
- Συγκρίνετε τον αριθμό σειράς στο κάτω μέρος του οργάνου με τον αριθμό σειράς στο δελτίο αποστολής.
 Αναφέρετε αμέσως οποιαδήποτε διαφορά.
- Ελέγξτε τα παρελκόμενα του οργάνου σε σχέση με αυτά που αναγράφονται στο δελτίο αποστολής.
- Αποθηκεύεστε τα υλικά συσκευασίας για τυχόν μεταφορά του οργάνου στο μέλλον.



3.2 Απαιτήσεις τροφοδοσίας

Το όργανο ανιχνεύει αυτόματα την παρεχόμενη τάση. Κατά συνέπεια, δεν είναι απαραίτητη οποιαδήποτε αλλαγή στο εύρος τάσης. Ελέγξτε τις προδιαγραφές τάσης και βεβαιωθείτε ότι η παρεχόμενη τάση είναι η σωστή σύμφωνα με τις παρακάτω προδιαγραφές:

Τάση:

-	
Βασικό όργανο με προσαρμογέα ΑC:	100 – 240 V AC, 50/60 Hz
Βασικό όργανο χωρίς προσαρμογέα ΑC:	24 V DC

Εάν η παραπάνω αναφερόμενη τάση δεν είναι διαθέσιμη στη χώρα σας, επικοινωνήστε με το τοπικό τμήμα εξυπηρέτησης πελατών της Tecan.

Συνδέετε το όργανο μόνο σε παροχή τροφοδοσίας με γείωση.





ΠΡΟΣΟΧΗ



ΠΡΟΣΟΧΗ

ΜΗΝ ΑΝΤΙΚΑΘΙΣΤΑΤΕ ΑΠΟΣΠΩΜΕΝΑ ΚΑΛΩΔΙΑ ΤΡΟΦΟΔΟΣΙΑΣ ΡΕΥΜΑΤΟΣ ΜΕ ΚΑΛΩΔΙΑ ΠΟΥ ΕΧΟΥΝ ΑΝΕΠΑΡΚΗ ΟΝΟΜΑΣΤΙΚΗ ΤΙΜΗ ΙΣΧΥΟΣ.

3.3 Περιβαλλοντικές απαιτήσεις

Το όργανο θα πρέπει να τοποθετείται σε επίπεδη επιφάνεια, στην οποία δεν υπάρχει σκόνη, διαλύτες και όξινοι ατμοί.

Για να εξασφαλίσετε σωστά αποτελέσματα, θα πρέπει να αποφεύγονται οι κραδασμοί και η άμεση έκθεση στο ηλιακό φως.

Θερμοκρασία περιβάλλοντος:	
Λειτουργία	15 °C έως 35 °C (59 °F έως 95 °F)
Αποθήκευση	-30 °C έως 60 °C (-22 °F έως 140 °F)
Σχετική υγρασία:	20% έως 80% χωρίς συμπύκνωση σε θερμοκρασία λειτουργίας



3.4 Απαιτήσεις συστήματος

	Υποστηρίζεται	Συνιστάται
Υπολογιστής	Υπολογιστής που να υποστηρίζει Windows με επεξεργαστή συμβατό με Pentium που τρέχει στο 1 GHz (διπύρηνος)	2 GHz (διπύρηνος)
Λειτουργικό	Windows 10 (32 bit)	
σύστημα	Windows 10 (64 bit)	
	Εκδόσεις: Ρro	
Μνήμη	Windows 10 (32-bit): 1 GB RAM	2 GB RAM
	Windows 10 (64-bit): 2 GB RAM	4 GB RAM
Ελεύθερος χώρος στο σκληρό δίσκο	3 GB	5 GB
Οθόνη	Κάρτα γραφικών Super VGA	
Ανάλυση	1024 x 600 και άνω	1920 x 1080
Βάθος χρώματος	256	
Ποντίκι	Ποντίκι της Microsoft ή συμβατή συσκευή κατάδειξης	
Επικοινωνία	1 x USB 2.0	2 x USB 2.0 1 x RS232 (σειριακή)
Συσκευές	Windows 10:	
	Συσκευή γραφικών DirectX 9 με WDDM 1.0 ή νεότερο πρόγραμμα οδήγησης	
.NET	Microsoft .NET Framework 3.5: Στα Windows 10 ο χρήστης θα κληθεί να εγκαταστήσει το απαιτούμενο πλαίσιο .NET (3.5), εάν δεν υπάρχει ήδη.	
Windows Installer	3.1	
	Εάν αυτή η έκδοση δεν υπάρχει, το πρόγραμμα εγκατάστασης/αναβάθμισης θα την εγκαταστήσει.	
Microsoft Excel	2007, 2010, 2013, 2016 (32 bit),	2010 (32 bit)
	2019 (32 bit),	2019 (32 bit)
	Microsoft Excel 365 (32 bit)	
	τποστηριζονται μονο εκδοσεις 32 bit!	
	ΔΕΝ υπου πριζονται εκοοσείς Starter!	

3.5 Ενεργοποίηση του οργάνου

Στις παρακάτω διαδικασίες περιγράφονται αναλυτικά τα απαραίτητα βήματα για την ενεργοποίηση του οργάνου.



ΠΡΟΣΟΧΗ

ΠΡΙΝ ΤΗΝ ΕΓΚΑΤΑΣΤΑΣΗ ΚΑΙ ΕΝΕΡΓΟΠΟΙΗΣΗ ΤΟΥ ΟΡΓΑΝΟΥ, ΑΥΤΟ ΘΑ ΠΡΕΠΕΙ ΝΑ ΕΧΕΙ ΤΟΠΟΘΕΤΗΘΕΙ ΟΡΘΙΟ ΤΟΥΛΑΧΙΣΤΟΝ ΓΙΑ ΤΡΕΙΣ ΩΡΕΣ ΩΣΤΕ ΝΑ ΑΠΟΦΕΥΧΘΕΙ Η ΣΥΜΠΥΚΝΩΣΗ ΠΟΥ ΘΑ ΜΠΟΡΟΥΣΕ ΝΑ ΠΡΟΚΑΛΕΣΕΙ ΒΡΑΧΥΚΥΚΛΩΜΑ.

Όταν οι παραπάνω αναφερόμενες απαιτήσεις πληρούνται, η εγκατάσταση πραγματοποιείται με την παρακάτω διαδικασία:

- 1. Συνδέστε το όργανο στον εξωτερικό υπολογιστή με το καλώδιο USB.
- Βεβαιωθείτε ότι ο διακόπτης κύριας τροφοδοσίας στην αριστερή πλευρά του οργάνου είναι στη θέση OFF (Απενεργοποίηση).
- Εισαγάγετε το καλώδιο τροφοδοσίας στην υποδοχή της κύριας τροφοδοσίας στην αριστερή πλευρά.
- Ενεργοποιήστε το όργανο χρησιμοποιώντας το διακόπτη της κύριας τροφοδοσίας στην αριστερή πλευρά.

Το όργανο είναι έτοιμο για τη μέτρηση μικροπλακών αφού εγκατασταθεί το λογισμικό.



Σημείωση Πριν ξεκινήσετε τις μετρήσεις βεβαιωθείτε ότι η θέση μικροπλάκας Α1 έχει τοποθετηθεί σωστά.

Η μέτρηση μικροπλακών μπορεί να γίνει μόνο χωρίς καλύμματα.

Κλείστε το κάλυμμα του φορέα μεταφοράς πλάκας πριν ξεκινήσετε μια μέτρηση ώστε να αποφύγετε την επίδραση του περιβάλλοντος φωτός στα αποτελέσματα.



Σημείωση Πάντα να βγάζετε τη μικροπλάκα από τη συσκευή ανάγνωσης αμέσως μετά την ολοκλήρωση της μέτρησης.



3.6 Λογισμικό

3.6.1 Εισαγωγή / Προοριζόμενη χρήση

Το λογισμικό ελέγχου του οργάνου και ανάλυσης δεδομένων **Magellan** παρέχεται με το όργανο.

To Magellan είναι ένα **λογισμό ελέγχου της συσκευής ανάγνωσης και** ανάλυσης δεδομένων γενικής χρήσης για την ανάλυση των δεδομένων που προέρχονται από ελέγχους μικροπλακών με τη χρήση συσκευών μέτρησης Tecan.

Το Magellan είναι διαθέσιμο σε δύο εκδόσεις:

- Magellan και
- Magellan Tracker

Το λογισμικό **Magellan** προορίζεται για προσδιορισμούς τελικού σημείου, κινητικής και πολλαπλής σήμανσης με το όργανο INFINITE F50 PLUS σύμφωνα με την προοριζόμενη χρήση. Ανατρέξτε στο κεφάλαιο 2.1 Προοριζόμενη χρήση.

Το **Magellan Tracker** προσφέρει όλη την απαραίτητη λειτουργικότητα ώστε να συμμορφώνεται ως προς τον κανονισμό 21, CFR τμήμα 11 του FDA, συμπληρωματικά της λειτουργικότητας του Magellan.



Σημείωση

Είναι σημαντικό να λάβετε υπόψη σας ότι η σωστή εγκατάσταση του οργάνου και του λογισμικού Magellan δεν εξασφαλίζουν τη συμμόρφωση με νομοθεσία και απαιτήσεις. Θα πρέπει να τεκμηριωθούν οι αντίστοιχες πολιτικές που αφορούν διεργασίες και τυπικές διαδικασίες λειτουργίας, συμπεριλαμβανομένων της επικύρωσης και του ελέγχου ποιότητας.

3.6.2 Διαδικασία εγκατάστασης

Για την εγκατάσταση του λογισμικού, τοποθετήστε το USB stick στη θύρα USB και ακολουθήστε τα παρακάτω βήματα:

- Ο οδηγός εγκατάστασης του Magellan θα πρέπει να ξεκινήσει αυτόματα και να σας καθοδηγήσει στη διαδικασία εγκατάστασης. Εάν δεν ξεκινήσει, πραγματοποιήστε εκτέλεση του αρχείου 'E:\Tecan.exe' (όπου Ε είναι το γράμμα της μονάδας οδήγησης του USB stick).
- Επιλέξτε 'Magellan software' (Λογισμικό Magellan) και ανάλογα με την έκδοση που έχετε προμηθευτεί - 'Install Magellan' (Εγκατάσταση Magellan) ή 'Install Magellan Tracker' (Εγκατάσταση Magellan Tracker) για να ξεκινήσετε τη διαδικασία εγκατάστασης. Κατόπιν, ακολουθήστε τον οδηγό εγκατάστασης.
- Κάντε κλικ στο Install (Εγκατάσταση) για να ξεκινήσετε τη διαδικασία εγκατάστασης του λογισμικού.
- Κάντε κλικ στο I accept the terms of the license agreement (Αποδέχομαι τους όρους του συμφωνητικού άδειας χρήσης) και στο Next (Επόμενο) για να συνεχίσετε.
- 5. Εμφανίζεται η σελίδα Customer Information (Πληροφορίες πελάτη): εισαγάγετε το όνομα χρήστη και τον οργανισμό.
- 6. Εμφανίζεται η σελίδα Configuration (Διαμόρφωση): επιλέξτε τη γλώσσα.
- Σελίδα Use for regulated environments (Χρήση σε ρυθμιζόμενα περιβάλλοντα): κάντε κλικ στο Next (Επόμενο) για να συνεχίσετε.



- 8. Κάντε κλικ στο Install (Εγκατάσταση) για να ξεκινήσετε την εγκατάσταση.
- Κάντε κλικ στο Finish (Τέλος) για να ολοκληρώσετε την εγκατάσταση και να κλείσετε το πρόγραμμα ρύθμισης.

Μπορείτε να ξεκινήσετε τη λειτουργία του λογισμικού από το μενού **Start** (Έναρξη) των Windows, επιλέγοντας **Magellan** στην ομάδα προγραμμάτων 'Tecan'.





Σημείωση Είναι πολύ σημαντικό ο άνθρωπος που θα πραγματοποιήσει την εγκατάσταση του λογισμικού να έχει δικαιώματα διαχειριστή στον υπολογιστή.

Σημείωση Η έκδοση V7.5 Magellan δεν μπορεί να εγκατασταθεί μαζί με άλλες εκδόσεις του λογισμικού Magellan.

Από προεπιλογή, όλοι οι τύποι αρχείων που σχετίζονται με το Magellan αποθηκεύονται στους αντίστοιχους υποκαταλόγους στον παρακάτω κατάλογο:

- 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 Επαλήθευση εγκατάστασης

Μπορείτε να ελέγξετε την επιτυχή εγκατάσταση του Magellan με το πρόγραμμα αυτόματης επαλήθευσης εγκατάστασης:

Πραγματοποιήστε εκτέλεση του αρχείου *TecanlQ.exe* από την προεπιλεγμένη διαδρομή εγκατάστασης (C:*Program Files\TecanWagellan) ή από το μενού* Start (Έναρξη) των Windows: Start (Έναρξη) > Programs (Προγράμματα) > Tecan > MagellanlQ.

Κάντε κλικ στο **Check** (Έλεγχος) για να ξεκινήσετε την επαλήθευση της εγκατάστασης. Όλα τα εγκατεστημένα στοιχεία θα πρέπει να έχουν ως κατάσταση το **OK**. Σε περίπτωση που αναφερθεί πρόβλημα, επικοινωνήστε με τον τοπικό αντιπρόσωπο πωλήσεων της Tecan.

Για να κλείσετε το πρόγραμμα επαλήθευσης της εγκατάστασης, κάντε κλικ στο **Cancel** (Άκυρο) ή **Exit** (Έξοδος).



Σημείωση Η επαλήθευση της εγκατάστασης θα πρέπει να επαναλαμβάνεται κάθε φορά που πραγματοποιείται εγκατάσταση ή ενημέρωση σε νεότερη έκδοση του Magellan.



3.6.4 Τα πρώτα βήματα με το Magellan

Ο βασικός τύπος περιβάλλοντος χρήσης στο Magellan είναι ο οδηγός.

Οι τυπικοί οδηγοί στο Magellan αντιπροσωπεύουν υποσυστήματα ροής εργασιών, τα οποία είναι «βήμα-προς-βήμα» οδηγοί για την εκτέλεση σύνθετων διαδικασιών.

Μερικές φορές υπάρχουν μενού, διαθέσιμα στη γραμμή κεφαλίδας. Τα Menu (Μενού) αποτελούν έναν κλασικό τρόπο χρήσης του λογισμικού:

το σχετικό στοιχείο μενού επιλέγεται από τα κύρια μενού.

Όλες οι επακόλουθες ενέργειες ξεκινούν άμεσα ή εμφανίζεται ένα πλαίσιο διαλόγου όταν υπάρχει η δυνατότητα πραγματοποίησης περαιτέρω επιλογών ή εισαγωγής στοιχείων.

Περιβάλλον χρήσης – Λίστα οδηγών

Μετά την εκκίνηση του Magellan, εμφανίζεται η οθόνη **Wizard List** (Λίστα οδηγών) :

Vizard List - Magellan		
\wedge	What do you want to do?	
Start measurement Evaluate result	 The Start Measurement wizard helps you to perform a measurement. You can either use a method or obtain raw data 	
Attach signature	e O	
Create/edit a method		
/		
Exit Magellan		

Η εκκίνηση κάθε οδηγού μπορεί να γίνει είτε με διπλό κλικ είτε επιλέγοντάς τον

		5
και κάνοντας κλικ στο κουμπί Next (Επόμενο)	Ľ	

Οδηγός έναρξης μέτρησης

Ο οδηγός Start Measurement (Έναρξη μέτρησης) περιλαμβάνει τις παρακάτω επιλογές:

- Η επιλογή Obtain Raw Data (Λήψη πρωτογενών δεδομένων) χρησιμοποιείται για τη γρήγορη και εύκολη δημιουργία πρωτογενών δεδομένων, με ρύθμιση των απαιτούμενων παραμέτρων μέτρησης και έναρξη μιας μέτρησης.
- Η επιλογή Use Predefined Method (Χρήση προκαθορισμένης μεθόδου) χρησιμοποιείται για την εκτέλεση μετρήσεων με βάση προκαθορισμένες μεθόδους.
- Η επιλογή Start Favorite (Έναρξη προτιμώμενης) χρησιμοποιείται για την επιλογή της πιο συχνά χρησιμοποιούμενης μεθόδου από τη λίστα με τα αριθμημένα εικονίδια.

Μετά την ολοκλήρωση της μέτρησης δημιουργείται ένα αρχείο χώρου εργασίας.



Οδηγός αξιολόγησης αποτελεσμάτων

Ο οδηγός **Evaluate Results** (Αξιολόγηση αποτελεσμάτων) χρησιμοποιείται για την προβολή των πρωτογενών δεδομένων και την αξιολόγηση των αποτελεσμάτων. Είναι δυνατή η προβολή των παραμέτρων αξιολόγησης και η επαναξιολόγηση των δεδομένων.

Οδηγός επισύναψης υπογραφής

Ο οδηγός **Attach Signature** (Επισύναψη υπογραφής) χρησιμοποιείται για την υπογραφή μεθόδων και αρχείων χώρου εργασίας. Αυτή η λειτουργία είναι διαθέσιμη μόνο στο Magellan Tracker.

Οδηγός δημιουργίας/επεξεργασίας λίστας αναγνωριστικών δειγμάτων

Ο οδηγός Create/Edit a Sample ID list (Δημιουργία/επεξεργασία λίστας αναγνωριστικών δειγμάτων) χρησιμοποιείται για τη δημιουργία νέων ή την επεξεργασία υπαρχουσών λιστών αναγνωριστικών δειγμάτων.

Οδηγός δημιουργίας/επεξεργασίας μεθόδου

Ο οδηγός **Create/Edit a method** (Δημιουργία/επεξεργασία μεθόδου) χρησιμοποιείται για τον ορισμό ή την επεξεργασία μεθόδων.



Σημείωση Για αναλυτικές πληροφορίες σχετικά με το λογισμικό, ανατρέξτε στις Οδηγίες Χρήσης Magellan.

Λάβετε υπόψη σας ότι μερικές λειτουργίες που περιγράφονται στις Οδηγίες Χρήσης Magellan ενδέχεται να μην είναι διαθέσιμες στην έκδοση V7.5 Magellan σε συνδυασμό με το INFINITE F50 PLUS. Ωστόσο, όλες οι απαραίτητες πληροφορίες περιγράφονται στον παρόν εγχειρίδιο.



Σημείωση Μπορείτε να βρείτε αναλυτικό παράδειγμα μέτρησης ELISA στο κεφάλαιο 6 Παράδειγμα εφαρμογής.

Αρχεία μεθόδου Sunrise που έχουν δημιουργηθεί με το Magellan V7.x ή νεότερη έκδοση μπορούν να ανοιχτούν με το Magellan V7.5. Οι παράμετροι μετρήσεων μετατρέπονται αυτόματα. Μέθοδοι Sunrise που έχουν δημιουργηθεί με προηγούμενες εκδόσεις πρέπει να μετατραπούν με τη χρήση της επιλογής 'Convert From' (Μετατροπή από) στο μενού miscellaneous/file handling (διάφορα/χειρισμός αρχείων).



3.7 Magellan – Πρόγραμμα επεξεργασίας παραμέτρων μέτρησης

Το Measurement Parameter Editor (Πρόγραμμα επεξεργασίας παραμέτρων μέτρησης) χρησιμοποιείται για τη ρύθμιση ροών εργασιών. Κάθε ροή εργασιών δημιουργείται εύκολα με τη μεταφορά και απόθεση των βημάτων της διεργασίας σε μια ακολουθία, σύμφωνα με την εφαρμογή. Κατόπιν, η ροή εργασιών της εφαρμογής είναι ορατή στο χρήστη από το παράθυρο ροής εργασιών. Κάθε βήμα της διεργασίας (στοιχείο του προγράμματος) μπορεί να αντιγραφεί και επικολληθεί (με τη χρήση των τυπικών συντομεύσεων των Windows Ctrl-C, Ctrl-V ή το σχετικό με το περιεχόμενο μενού του ποντικιού) και να μετακινηθεί στην επιθυμητή θέση μέσα στη ροή εργασιών.

Measurement parameters - infinite F5	0		
🔶 Lab Ware 🔹	🐟 ▼ Plate	1	Selection
Interview Plate	Dian definition (ICDE 001) - Conient 02 Ella Transmand	Detailt	Nothing selected
🚸 Part of Plate	nae deimion. [dinc bok] - diemei bo nati nanspalenk	Use a part of the plate	
Measurements	A D - 2011		
Absorbance	PratorPlate ALHI2		
	Absorbance	3	
Actions (8)	Wavelength Measurement 405 mm		
Nove Plate	Reference: 405 nm		
Kinetic S			
A Kinetic Condition			
Priscellaneous			
 User Request 			
Wait (Timer)			
J Incubation			
			Παράθυρο
Γοαμμή ελένχου	νώιτρηνος ρομθάρη		πληροφο-
. beinheit enertVee			
			ριών
		~	
Number of plates: 1			
Help			_
Cancel (<<< Back	CHOOSE M	EASUREMENT PARAMET	ERS
© 2009 Tecan			

Το **Measurement Parameter Editor** (Πρόγραμμα επεξεργασίας παραμέτρων μέτρησης) αποτελείται από τα παρακάτω στοιχεία, τα οποία περιγράφονται αναλυτικά στα επόμενα κεφάλαια:

- Γραμμή ελέγχου
- Παράθυρο ροής εργασιών
- Παράθυρο πληροφοριών



3.7.1 Γραμμή ελέγχου

Η **Γραμμή ελέγχου** χωρίζεται σε πέντε τμήματα. Κάθε τμήμα περιέχει στοιχεία προγράμματος που χρησιμοποιούνται για τη δημιουργία μιας μεμονωμένης ροής εργασιών.

Μπορείτε να δημιουργήσετε μια ροή εργασιών είτε με διπλό κλικ στο επιλεγμένο στοιχείο προγράμματος είτε με μεταφορά και απόθεσή του στο παράθυρο ροής εργασιών.

Κατά τη χρήση του INFINITE F50 PLUS είναι διαθέσιμα τα παρακάτω στοιχεία προγράμματος:

Lab Ware (Εργαστήριο)	Plate (Πλάκα) Part of Plate (Τμήμα πλάκας)	
Measurements (Μετρήσεις)	Absorbance (Απορρόφηση)	
Actions (Ενέργειες) Shaking (Ανακίνηση) Μονe Plate (Μετακίνηση πλάκας)		
Kinetic (Ανάλυση κινητικής)	νάλυση Kinetic Cycle (Κύκλος ανάλυσης κινητικής) Kinetic Condition (Συνθήκη ανάλυσης κινητικής)	
Miscellaneous (Διάφορα)	Comment (Σχόλιο) User Request (Αίτημα χρήστη) Wait (Timer) (Αναμονή [Χρονοδιακόπτης]) Incubation (Επώαση)	

Εργαστήριο

Πλάκα

Το στοιχείο προγράμματος **Plate** (Πλάκα) χρησιμοποιείται για την επιλογή μορφής πλάκας από την αναπτυσσόμενη λίστα **Plate definition** (Ορισμός πλάκας). Κάντε κλικ στο **Details**...(Λεπτομέρειες) για να δείτε περισσότερες πληροφορίες για την επιλεγμένη πλάκα.

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

Τμήμα πλάκας

Από προεπιλογή, το στοιχείο προγράμματος **Part of Plate** (Τμήμα πλάκας)είναι συμπτυγμένο. Κάνοντας κλικ στο αναπτύσσεται και εμφανίζει μια μικροπλάκα 96 κυψελίδων. Για να μετρήσετε μεμονωμένες κυψελίδες, κάντε κλικ στην κυψελίδα που θέλετε. Για να μετρήσετε ένα εύρος κυψελίδων, σύρετε ένα πλαίσιο γύρω από το επιθυμητό εύρος. Κάνοντας κλικ στο **Details...** (Λεπτομέρειες), μπορείτε να κάνετε ζουμ στην προεπισκόπηση της πλάκας.

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



Ανεξάρτητα τμήματα πλάκας

Μπορείτε να επιλέξετε ανεξάρτητα τμήματα πλάκας:



Μπορείτε να επιλέξετε ένα δεύτερο εύρος κυψελίδων πατώντας το πλήκτρο Control στο πληκτρολόγιο και σύροντας ένα πλαίσιο πάνω από τις προς επιλογή κυψελίδες.

Μετρήσεις

Απορρόφηση

Το στοιχείο προγράμματος **Absorbance** (Απορρόφηση) χρησιμοποιείται για την πραγματοποίηση μετρήσεων απορρόφησης. Εισαγάγετε ή επιλέξτε τις αντίστοιχες παραμέτρους.

Τα διαθέσιμα μήκη κύματος μετρήσεων και φίλτρων αναφοράς εμφανίζονται σε δύο αναπτυσσόμενες λίστες, σύμφωνα με τα φίλτρα απορρόφησης που έχουν τοποθετηθεί. Εάν οι αναπτυσσόμενες λίστες είναι άδειες, τα φίλτρα δεν έχουν οριστεί.

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

Ενέργειες

Ανακίνηση

Επιλέξτε το στοιχείο προγράμματος **Shaking** (Ανακίνηση) εάν η πλάκα πρέπει να ανακινηθεί, είτε πριν τη μέτρηση είτε μεταξύ των κύκλων ανάλυσης κινητικής.

📉 🔻 Shaking		4
Parameter Duration: 1 📚 sec Intensity: V A	Wide vmplitude: 14,1 mm irequency: 2,1 Hz	Wait a couple of seconds



3. Τα πρώτα βήματα

Εισαγάγετε τις αντίστοιχες παραμέτρους:

Duration (Διάρκεια)	Εισαγάγετε τη διάρκεια της διαδικασίας ανακίνησης.
Intensity (Ένταση)	Εισαγάγετε την επιθυμητή λειτουργία ανακίνησης. Κατά την επιλογή της αντίστοιχης λειτουργίας ανακίνησης, εμφανίζονται το εύρος και η συχνότητα.

Για τις διαθέσιμες λειτουργίες ανακίνησης, ανατρέξτε στο κεφάλαιο 4.1.1 Ανακίνηση μικροπλάκας.

Κάνοντας κλικ στη σύνδεση <u>Wait a couple of seconds</u> (Αναμονή λίγα δευτερόλεπτα) μπορείτε να εισάγετε ένα νέο στοιχείο προγράμματος. Για λεπτομέρειες, ανατρέξτε στη σελίδα 25.

Ανάλυση κινητικής

Κύκλος ανάλυσης κινητικής

Χρησιμοποιήστε το στοιχείο προγράμματος **Kinetic Cycle** (Κύκλος ανάλυσης κινητικής)**για την πραγματοποίηση πολλαπλών συνεχόμενων μετρήσεων**, οι οποίες μπορεί να πραγματοποιηθούν σε συγκεκριμένα χρονικά διαστήματα.

🎧 🔻 Kinetic Cycle		3
Oycles Image: Number of cycles: Duration	Kinetic Interval ✓ Use kinetic interval: ● Time: 00:01:00 ♀ (hh:mm:ss) ● Time: 60000 ♦	

Εισαγάγετε τις αντίστοιχες παραμέτρους:

Cycles (Κύκλοι)	Number of cycles (Αριθμός κύκλων): Εισαγάγετε έναν αριθμό ή κάντε κλικ στα βελάκια πάνω ή κάτω για τον αριθμό των βημάτων μέτρησης (2 - 1000 κύκλοι) Duration (Διάρκεια): Εισαγάγετε την διάρκεια με τη μορφή ωω:λλ:δδ.
Kinetic Interval	Use kinetic interval (Χρήση χρονικού διαστήματος
(Χρονικό διάστημα	ανάλυσης κινητικής): Εισαγάγετε το χρονικό διάστημα
ανάλυσης κινητικής)	(ωω:λλ:δδ ή ms).

Συνθήκη ανάλυσης κινητικής

Χρησιμοποιήστε το στοιχείο προγράμματος **Kinetic Condition** (Συνθήκη ανάλυσης κινητικής) για να ορίσετε ποιες ενέργειες θα πραγματοποιηθούν σε συγκεκριμένο κύκλο.

🙊 🔻 Kinetic Condition	4
Execute commands at cycle:	

Εάν **στο Execute command at cycle** (Εκτέλεση εντολής στον κύκλο) εισάγετε το **2**, εντός μιας μέτρησης ανάλυσης κινητικής, που περιέχει π.χ. ένα βήμα **Shake** (Ανακίνηση), η ανακίνηση πραγματοποιείται μόνο στον κύκλο 2.



(\mathbf{i})

3. Τα πρώτα βήματα

Σημείωση

Οι συνθήκες ανάλυσης κινητικής, όπως η ανακίνηση, θα πρέπει να εισαχθούν αμέσως μετά από ένα στοιχείο προγράμματος Kinetic Cycle (Κύκλος κινητικής ανάλυσης) ώστε να εξασφαλιστεί η βέλτιστη δυνατότητα αναπαραγωγής των αποτελεσμάτων. Προτείνεται στους χρήστες η δημιουργία κατάλληλων δεσμών ενεργειών πριν από τις μετρήσεις και η χρήση της ίδιας δέσμης για όλες τις παρόμοιες μετρήσεις ανάλυσης κινητικής με στόχο τη λήψη συγκρίσιμων αποτελεσμάτων.

Διάφορα

Σχόλιο

Χρησιμοποιήστε το στοιχείο προγράμματος **Comment** (Σχόλιο) για να εισάγετε στο πεδίο κείμενου ένα σχόλιο ή παρατήρηση για την τρέχουσα μέτρηση.

💷 🔻 Comment	6
Courses I	

Αίτημα χρήστη

Το στοιχείο προγράμματος **User Request (Αίτημα χρήστη)** πληροφορεί το χειριστή του οργάνου να εκτελέσει μια συγκεκριμένη ενέργεια κατά τη ροή εργασιών, σε συγκεκριμένη στιγμή.

🔷 🔻 User Request	7
Text	

Εάν, για παράδειγμα, το στοιχείο προγράμματος **Move Plate** (Μετακίνηση πλάκας) χρησιμοποιείται για τη μετακίνηση της πλάκας εκτός, για την εκτέλεση μιας συγκεκριμένης ενέργειας, το κείμενο που θα εισαχθεί θα πρέπει να πληροφορεί το χειριστή να εκτελέσει αυτές τις ενέργειες. Ένα πλαίσιο διαλόγου εμφανίζει το μήνυμα και η διαδικασία μέτρησης διακόπτεται έως ότου κάνετε κλικ στο **OK**.

Αναμονή (Χρονοδιακόπτης)

Χρησιμοποιήστε το στοιχείο προγράμματος **Wait (Timer)** (Αναμονή [Χρονοδιακόπτης]) για να ορίσετε μια συγκεκριμένη περίοδο αναμονής πριν την εκτέλεση του επόμενου βήματος σε μια ροή εργασιών.

Στο πεδίο Wait time (Χρόνος αναμονής), εισαγάγετε τον απαιτούμενο χρόνο.

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



3. Τα πρώτα βήματα

Εισαγάγετε τις αντίστοιχες παραμέτρους:

Timer (Χρονοδιακόπτης)	Εισαγάγετε Wait time (Χρόνος αναμονής) (ωω:λλ:δδ)
Options (Επιλογές)	Ignore wait at last kinetic cycle (Να αγνοηθεί η αναμονή κατά τον τελευταίο κύκλο ανάλυσης κινητικής): Όταν το βήμα προγράμματος Wait (Timer) (Αναμονή [Χρονοδιακόπτης]) είναι η τελευταία ενέργεια σε έναν κύκλο ανάλυσης κινητικής, ο χρόνος αναμονής θα αγνοηθεί στον τελευταίο κύκλο.

Επώαση

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

Εισαγάγετε τις κατάλληλες παραμέτρους για την επώαση:

Incubation time	Εισαγάγετε το συνολικό χρόνο (τουλάχιστον 5
(Χρόνος επώασης)	δευτερόλεπτα)

3.7.2 Παράθυρο ροής εργασιών

Το Παράθυρο ροής εργασιών είναι το παράθυρο, όπου είναι ορατή η δέσμη ενεργειών μέτρησης και όπου πραγματοποιείται ο ορισμός και η επεξεργασία των παραμέτρων.

Υπάρχουν δύο τρόποι με τους οποίους μπορείτε να εισάγετε ένα στοιχείο προγράμματος από τη **Γραμμή ελέγχου** στο **Παράθυρο ροής εργασιών**:

- Επιλέξτε ένα στοιχείο προγράμματος από τη Γραμμή ελέγχου. Κάνοντας διπλό κλικ πάνω σε αυτό πραγματοποιείται η εισαγωγή του στο Παράθυρο ροής εργασιών αμέσως μετά το προηγούμενο στοιχείο προγράμματος.
- Κάντε κλικ στο στοιχείο προγράμματος στη Γραμμή ελέγχου και σύρετέ το στο Παράθυρο ροής εργασιών στην αντίστοιχη θέση.

Τα στοιχεία προγράμματος αριθμοδοτούνται σύμφωνα με την ακολουθία τους.

Μετά την εισαγωγή ενός στοιχείου προγράμματος στο Παράθυρο ροής εργασιών, μπορείτε να εισάγετε ή επεξεργαστείτε τις ρυθμίσεις και τις παραμέτρους αυτού του στοιχείου.

Τα στοιχεία προγράμματος μέσα στο Παράθυρο ροής εργασιών μπορούν να συμπτυχθούν ώστε να εμφανίζονται οι πιο σημαντικές πληροφορίες ή να αναπτυχθούν ώστε να υπάρχει πρόσβαση σε όλες τις επεξεργάσιμες λειτουργίες. Κάντε κλικ σε ένα από τα τρίγωνα δίπλα στον τίτλο ενός στοιχείου προγράμματος, ή , για εναλλαγή των δύο προβολών.

Από προεπιλογή, το πρόγραμμα επεξεργασίας παραμέτρων μέτρησης ξεκινά με το στοιχείο Plate (Πλάκα), το στοιχείο Part of Plate (Τμήμα πλάκας) (σε σύμπτυξη) και ένα στοιχείο Absorbance (Απορρόφηση) στο Παράθυρο ροής εργασιών.

Τα τρέχοντα επιλεγμένα στοιχεία προγράμματος μέσα στο Παράθυρο ροής εργασιών εμφανίζονται με μια κίτρινη γραμμή στην πάνω πλευρά.

Εάν ένα στοιχείο προγράμματος περιέχει σφάλματα ή δεν είναι έγκυρο μέσα στην τρέχουσα ροή εργασιών, το στοιχείο θα σημανθεί με ένα σημάδι σφάλματος και ο αριθμός του στοιχείου επισημαίνεται με κόκκινο. Στο Παράθυρο πληροφοριών, εμφανίζονται αναλυτικές πληροφορίες για το σφάλμα. Εάν η ροή εργασιών περιέχει σφάλματα, δεν είναι δυνατή η επιλογή παραμέτρων μέτρησης.



Ιεραρχία στοιχείων

Η ιεραρχία των στοιχείων στο Παράθυρο ροής εργασιών έχει ως εξής:

- 1. Πλάκα
- 2. Τμήμα πλάκας (Εύρος)

Όποιο βήμα μέτρησης θέλετε μπορεί να εισαχθεί αμέσως μετά από ένα στοιχείο πλάκας ή εύρους. Χρησιμοποιήστε τις επιλογές **Release** (Αποδέσμευση) και **Indent** (Δέσμευση) για να τροποποιήσετε την ακολουθία εκτέλεσης ενός τμήματος λωρίδας. Επιλέξτε ένα στοιχείο στο **Παράθυρο ροής εργασιών**, κάντε δεξί κλικ και επιλέξτε **Release** (Αποδέσμευση) ή **Indent** (Δέσμευση).

Μπορείτε να εισάγετε άλλα στοιχεία από τη **Γραμμή ελέγχου** στην ιεραρχία μιας ροής εργασιών, ως εξής:

Το πρώτο στοιχείο **Range** (Εύρος) μπορεί να εισαχθεί αμέσως μετά το στοιχείο **Plate** (Πλάκα). Κατόπιν, μπορείτε να εισάγετε όλα τα επόμενα στοιχεία **Range** (Εύρος).

Είναι επίσης δυνατή η χρήση βημάτων ανάλυσης κινητικής μέσα σε ένα στοιχείο **Plate** (Πλάκα) ή **Range** (Εύρος).

Τα βήματα User Request (Αίτημα χρήστη), **Comment** (Σχόλιο) και **Wait** (Αναμονή) είναι εφικτά μέσα σε ένα στοιχείο **Plate** (Πλάκα) ή **Range** (Εύρος).

3.7.3 Παράθυρο πληροφοριών

Το Παράθυρο πληροφοριών στη δεξιά πλευρά της οθόνης εμφανίζει πληροφορίες που σχετίζονται με το τρέχον επιλεγμένο στοιχείο προγράμματος. Τυχόν προειδοποιήσεις ή σφάλματα εμφανίζονται εδώ.

3.8 Magellan – Ορισμός μετρήσεων

Στο παρακάτω κεφάλαιο περιγράφονται μερικά παραδείγματα για τον ορισμό διαφορετικών μετρήσεων.

3.8.1 Ορισμός μετρήσεων τελικού σημείου

Το παρακάτω παράδειγμα περιγράφει τη λειτουργία **Absorbance Endpoint Measurement** (Μέτρηση τελικού σημείου απορρόφησης)σε όλες τις κυψελίδες μιας μικροπλάκας 96 κυψελίδων.

- Επιλέξτε μια μικροπλάκα 96 κυψελίδων από την αναπτυσσόμενη λίστα Plate definition (Ορισμός πλάκας).
- Από προεπιλογή, όλες οι κυψελίδες της μικροπλάκας 96 κυψελίδων είναι επιλεγμένες προς μέτρηση.
- 3. Εισαγάγετε τα επιθυμητά μήκη κύματος μέτρησης και αναφοράς.

Part of Plate 1 2 3 4 5 6 7 8 9 10 11 12 A B C	I 2 3 4 5 6 7 8 9 10 11 12 I 2 3 4 5 6 7 8 9 10 11 12 Details I 0	Plate definition:	[COS96it] - Coming 96 Flat Transparent	Use a part of the pla
✓ Absorbance Wavelength Messurement: 450 nm ✓ Reference: 620 nm	Vavelength Messurement: 450 nm Varme Label Reference: 620 nm V		Part of Plate	
		M		
			Reference: 620 nm	
	/			~



3.8.2 Ορισμός μετρήσεων πολλαπλής σήμανσης

Το παρακάτω παράδειγμα περιγράφει μια μέτρηση **Absorbance Multilabel Measurement** (Μέτρηση απορρόφησης πολλαπλής σήμανσης)σε ένα καθορισμένο εύρος από μικροπλάκες 96 κυψελίδων (A1:E7). Θα μετρηθούν τρεις σημάνσεις απορρόφησης.

- Επιλέξτε μια μικροπλάκα 96 κυψελίδων από την αναπτυσσόμενη λίστα Plate definition (Ορισμός πλάκας).
- Από προεπιλογή, όλα τα φρεάτια της μικροπλάκας 96 κυψελίδων είναι επιλεγμένα προς μέτρηση.
 Κάντε κλικ στο για να αναπτύξετε το στοιχείο Part of Plate (Τμήμα πλάκας). Κατόπιν, επιλέξτε το επιθυμητό εύρος πλακών (A1:E7).
- 3. Εισαγάγετε το επιθυμητό μήκος κύματος μέτρησης.
- 4. Εισαγάγετε 2 ακόμα στοιχεία **Absorbance** (Απορρόφηση) και τα μήκη κύματος μέτρησης.

🔶 🔻 Plat	3	1
Plate definition	[COS96ft] - Corning 96 Flat Transparent	V Details
		Use a part of the plate
- 🚸 🔻	Part of Plate	2
A B C C E G G H	2 3 4 5 8 7 8 9 10 11 12 Detaits	
	▼ Absorbance	3
ſ	Wavelength	
	Measurement: 405 nm Vame: Label Reference: 405 nm V	
	▼ Absorbance	4
ſ	Wavelength Label	
	Measurement: 492 nm 💌 Name: Label	
	Reference: 405 nm	
	▼ Absorbance	5
	Wavelength Label	
	Measurement: 620 nm Vame: Label	
	Reference: 405 nm	
		- C



3.8.3 Ορισμός μετρήσεων ανάλυσης κινητικής

Το παρακάτω παράδειγμα περιγράφει μια μέτρηση ανάλυσης κινητικής μιας μικροπλάκας 96 κυψελίδων.

- Επιλέξτε μια μικροπλάκα 96 κυψελίδων από την αναπτυσσόμενη λίστα Plate definition (Ορισμός πλάκας).
- Εισαγάγετε ένα στοιχείο προγράμματος Kinetic Cycle (Κύκλος ανάλυσης κινητικής) μεταξύ του τμήματος της πλάκας και του στοιχείου απορρόφησης.
- 3. Cycles/Number of cycles (Κύκλοι/Αριθμός κύκλων): 50
- 4. Kinetic interval (Χρονικό διάστημα ανάλυσης κινητικής) (το χρονικό διάστημα μεταξύ μετρήσεων): επιλέξτε Use kinetic interval (Χρήση χρονικού διαστήματος ανάλυσης κινητικής) και εισαγάγετε: 2 λεπτά 30 δευτερόλεπτα.
- 5. Ορίστε το στοιχείο **Absorbance** (Απορρόφηση) εισάγοντας το επιθυμητό μήκος κύματος μέτρησης.

			Use a part of the plat
🔹 🔹 🔹	Part of Plate		:
A B C D C C C C C C C C C C C C C C C C C		etails	
G	🔻 Kinetic Cycle		:
-0	ycles	Kinetic Interval	
0	Duration	✓ Use kinetic interval: ⊙ Time: 00:02:30 ♀ ○ Time: 150000 ♀ ms	
	🖕 🔻 Absorbance		
	Wavelength	Label	
	Measurement: 492 nm	Name: Label1 🗸	
			~



3.8.4 Δέσμευση και αποδέσμευση στοιχείων προγράμματος

Η δέσμευση/αποδέσμευση ενός στοιχείου προγράμματος τροποποιεί τη ροή εργασιών του οργάνου κατά τη διάρκεια μετρήσεων.

Οι ενέργειες όλων των στοιχείων προγράμματος με τον ίδιο βαθμό δέσμευσης εκτελούνται διαδοχικά. Η μόνη εξάρτηση μεταξύ αυτών των στοιχείων προγράμματος είναι ότι η επόμενη ενέργεια ξεκινά αμέσως μόλις ολοκληρωθεί η προηγούμενη.

Ένα στοιχείο προγράμματος το οποίο έχει μεγαλύτερο βαθμό δέσμευσης από ό,τι το προηγούμενο στοιχείο προγράμματος υποδεικνύει εξάρτηση μεταξύ αυτών των δύο στοιχείων. Αυτό σημαίνει ότι οι παράμετροι που έχουν οριστεί στο πρώτο στοιχείο προγράμματος είναι επίσης ενεργές και για το δεύτερο (δεσμευμένο) στοιχείο.

Ακολουθεί ένα παράδειγμα για τον τρόπο ορισμού **ανάλυσης κινητικής** πολλαπλής σήμανσης με δύο σημάνσεις απορρόφησης. Αυτό το παράδειγμα δείχνει ότι τα δύο στοιχεία προγράμματος Absorbance (Απορρόφηση) εξαρτώνται από το στοιχείο προγράμματος Kinetic Cycle (Κύκλος κινητικής ανάλυσης), το οποίο εξαρτάται από το στοιχείο προγράμματος Part of Plate (Τμήμα πλάκας), το οποίο, με τη σειρά του, εξαρτάται από το στοιχείο προγράμματος Plate (Πλάκα). Ορίστε τις παραμέτρους για ένα παράδειγμα, ως εξής:

- 1. Plate (Πλάκα): π.χ. Greiner 96, επίπεδη, διαφανής
- Kinetic Cycle/Number of cycles (Κύκλος ανάλυσης κινητικής/Αριθμός κύκλων):
 5
- 3. Absorbance/Wavelength Label 1 (Απορρόφηση/Σήμανση μήκους κύματος 1): 450 nm
- Absorbance/Wavelength Label 2 (Απορρόφηση/Σήμανση μήκους κύματος 2): 492 nm



		_
Plate definition:	[GRE96it] - Greiner 96 Flat Transparent	Vetails
		Use a part of the pla
🔶 🛨	Part of Plate	
A B C C C C C C C C C C C C C C C C C C	2 3 4 5 6 7 8 9 10 11 12 Details Details	
G) ▼ Kinetic Cycle	
C.	Cycles Kinetic Interval	
	Vumber of cycles: 5 Duration Kinetic Interval	
	Ouration Kinetic Interval Duration Use kinetic interval Use kinetic interval Use kinetic interval	_
	Vamber of cycles: 5 Duration Use kinetic interval Variation Use kinetic interval Wavelength Label	
	Ouration Kinetic Interval Duration Use kinetic interval Wavelength Label Measurement: 450 nm Reference: 405 nm	
	Vamber of cycles: 5 Duration Use kinetic interval Use kinetic interval Wavelength Messurement: 450 nm Reference: 405 nm Wavelength Label Name: Label1 Vamber	
	Vamber of cycles: 5 Duration Use kinetic interval Wavelength Label Messurement: 450 nm Reference: 405 nm Wavelength Label Wavelength Label Wavelength Label	
	Sycles Number of cycles: 5 Duration Use kinetic interval Wavelength Label Messurement: 405 mm 405 mm Label Wavelength Label Wavelength Label Wavelength Label Wavelength Name: Label Varelength	
	Varber Number of cycles: 5 Duration Use kinetic interval Use kinetic interval Wavelength Reference: 405 nm Wavelength Name: Label Name: Label	

Το Παράθυρο ροής εργασιών εμφανίζεται όπως στην παρακάτω εικόνα:

Ο παραπάνω ορισμός έχει ως αποτέλεσμα την παρακάτω ροή εργασιών:

Η απορρόφηση όλων των κυψελίδων της μικροπλάκας των 96 κυψελίδων μετριέται πρώτα στα 450 nm και κατόπιν στα 492 nm. Και οι δύο μετρήσεις απορρόφησης πραγματοποιούνται σε 5 κύκλους ανάλυσης κινητικής.



Η αποδέσμευση του δεύτερου στοιχείου προγράμματος **Absorbance** (Απορρόφηση), έτσι ώστε να είναι ευθυγραμμισμένο με το στοιχείο **Kinetic Cycle** (Κύκλος ανάλυσης κινητικής), τροποποιεί τη ροή εργασιών. Επιλέξτε το δεύτερο στοιχείο προγράμματος **Absorbance** (Απορρόφηση) και κάντε δεξί κλικ. Επιλέξτε **Release Strip** (Αποδέσμευση λωρίδας) από το σχετικό με το περιεχόμενο μενού. Το **Παράθυρο ροής εργασιών** εμφανίζεται όπως στην παρακάτω εικόνα:

Plate definiti	ion: [GRE96it] - Greiner 96 Flat Transparent	V Details
		Use a part of the plat
-	▼ Part of Plate	:
A B C D F G H	1 2 3 4 5 6 7 8 9 10 11 12 0 <th></th>	
	😡 ▼ Kinetic Cycle	:
	Cycles Kinetic Interval	
	Opcles Signification O Duration Use kinetic interval	
	Cycles Signification Image: Signification of cycles: Signification Duration Use kinetic interval Image: Signification of cycles: Signification Image: Signification Signification </td <td></td>	
	Opcles 5 O Number of cycles: 5 Duration Use kinetic interval Image: the structure of cycles: 1 Image: the structure of cycles: 5	
	Cycles 5 • Number of cycles: 5 • Duration □ Use kinetic interval • Mayelength • Label • Measurement: • 450 nm • Label	_
	Cycles S • Number of cycles: Duration • Duration • Use kinetic interval • Wavelength • Label Measurement: 450 nm • Reference: • Optime	
	Cycles 5 Image: Solution Use kinetic interval Image: Duration Image: Use kinetic interval Image: Wavelength Image: Use kinetic interval Image: Reference: 405 nm Image: Use kinetic interval Image: Wavelength Image: Use kinetic interval<	
	Cycles Signification Image: Signification of cycles: Signification Duration Image: Use kinetic interval Image: Wavelength Label Measurement: 450 nm Image: Reference: 405 nm	
	Cycles Signification Image: Signification of cycles: Signification Image: Duration Image: Signification of cycles: Image: Maxelength Image: Label Image	
	Cycles Signification Image: Signification of cycles: Signification Duration Image: Signification of cycles: Image: Duration Image: Duration	
	Cycles Kinetic Interval Image: Signal Action of Cycles: Signal Action of Cycles: Image: Duration Image: Duration	

Σε αυτήν τη ροή εργασιών, μια μέτρηση **απορρόφησης ανάλυσης κινητικής** με 5 κύκλους εκτελείται για την πρώτη απορρόφηση στα 450 nm. Μετά την ολοκλήρωση αυτής της ακολουθίας, εκτελείται η μέτρηση **απορρόφησης τελικού σημείου** στα 492 nm.

3.9 Βελτιστοποίηση για καλύτερη απόδοση

Το όργανο έχει δοκιμαστεί πλήρως στο εργοστάσιο ώστε να εξασφαλίζεται ότι η απόδοσή του θα είναι εντός των προσδιορισμένων ορίων (για λεπτομέρειες, ανατρέξτε στο κεφάλαιο 4.4.2 Προδιαγραφές μέτρησης).

Η τήρηση των παρακάτω συστάσεων θα έχει ως αποτέλεσμα τη λειτουργία του οργάνου με τη μεγαλύτερη δυνατή ορθότητα.

3.9.1 Θέση οργάνου

Το όργανο θα πρέπει να τοποθετηθεί στην κατάλληλη θέση (για αναλυτικές πληροφορίες, ανατρέξτε στο κεφάλαιο 3.3 Ασφάλεια).



3.9.2 Διαδικασία λειτουργίας

Γενικά

- Συνιστάται η τήρηση των τυπικών διαδικασιών λειτουργίας για τους προσδιορισμούς που χρησιμοποιούνται.
- Η καλύτερη δυνατότητα αναπαραγωγής λαμβάνεται όταν το μήκος κύματος της μέτρησης αντιστοιχεί στο μέγιστο μήκος κύματος απορρόφησης του συγκεκριμένου διαλύματος.
 Είναι σημαντικό να χρησιμοποιείτε το μέγιστο μήκος κύματος απορρόφησης εάν η καμπύλη απορρόφησης του δείγματος είναι πάνω από ένα στενό φάσμα μήκους κύματος.
 Λάβετε υπόψη ότι οι μετρήσεις στην κλίση μιας κορυφής απορρόφησης θα περιορίσουν την ορθότητα των τιμών OD.
- Μετά την πραγματοποίηση μέτρησης κάθε πλάκας, ανατρέξτε στο πακέτο του κιτ ελέγχου για πληροφορίες σχετικά με τη διαδικασία επικύρωσης.
- Χρησιμοποιείτε τα συνιστώμενα φίλτρα απορρόφησης για το INFINITE F50 PLUS.

Μικροπλάκες

 Το όργανο μπορεί να χρησιμοποιηθεί με τους τύπους μικροπλακών που περιγράφονται στο κεφάλαιο 4.4.3 Μικροπλάκες. Τα καλύτερα αποτελέσματα λαμβάνονται όταν χρησιμοποιούνται μικροπλάκες επίπεδου πυθμένα. Ανάλογα με τον τύπο μικροπλάκας που χρησιμοποιείται, τα αποτελέσματα της μέτρησης ενδέχεται να διαφέρουν.

Να είστε ιδιαίτερα προσεκτικοί κατά τη χρήση μικροπλακών με σχήμα πυθμένα C, U ή V ή πλακών λωρίδων κυψελίδων καθώς είναι πιθανόν τα αποτελέσματα της μέτρησης να διαφέρουν σε σχέση με τις προδιαγραφές που περιγράφονται στο παρόν έγγραφο. Βεβαιωθείτε ότι ο τύπος μικροπλάκας που χρησιμοποιείται με το μετρητή απορρόφησης INFINITE F50 PLUS είναι κατάλληλος για την αντίστοιχη εφαρμογή.

- Χρησιμοποιείτε μόνο μικροπλάκες που είναι απόλυτα καθαρές.
- Αποφεύγετε την επικάθιση σκόνης πάνω στα διαλύματα ή τη μικροπλάκα κατά την περίοδο επώασης πριν την μέτρηση.
 Συνιστάται η χρήση καλύμματος για προστασία κάθε φορά που μια μικροπλάκα τοποθετείται εκτός του οργάνου.
- Οι ανακρίβειες στην ποσότητα διαλύματος που μεταφέρεται με πιπέτα επηρεάζουν περισσότερο τα αποτελέσματα που θα ληφθούν, όταν χρησιμοποιούνται μικρές ποσότητες διαλύματος.
- Η μορφή του μηνίσκου του διαλύματος μπορεί να προκαλέσει ανακρίβειες στα αποτελέσματα, ειδικά εάν χρησιμοποιούνται μικρές ποσότητες διαλύματος.



4. Λειτουργίες οργάνου

4.1 Λειτουργίες οργάνου

Οι παρακάτω λειτουργίες μέτρησης απορρόφησης είναι διαθέσιμες στο INFINITE F50 PLUS:

μετρήσεις τελικού σημείου, ανάλυσης κινητικής και πολλαπλής σήμανσης.

4.1.1 Ανακίνηση μικροπλάκας

To INFINITE F50 PLUS έχει τη δυνατότητα να ανακινήσει τη μικροπλάκα πριν τη μέτρηση. Επίσης, η μικροπλάκα μπορεί να ανακινηθεί μεταξύ των κύκλων μέτρησης ανάλυσης κινητικής.

Χρησιμοποιήστε το Magellan για τη ρύθμιση της λειτουργίας ανακίνησης.



ΠΡΟΣΟΧΗ ΚΑΤΑ ΤΗΝ ΑΝΑΚΙΝΗΣΗ ΜΙΚΡΟΠΛΑΚΩΝ, ΕΝΔΕΧΕΤΑΙ ΝΑ ΧΥΘΕΙ ΠΕΡΙΕΧΟΜΕΝΟ ΕΑΝ ΟΙ ΚΥΨΕΛΙΔΕΣ ΓΕΜΙΣΟΥΝ ΜΕ ΠΕΡΙΣΣΟΤΕΡΟ ΟΓΚΟ ΑΠΟ ΤΟ ΚΑΝΟΝΙΚΟ.

Λειτουργίες ανακίνησης για το INFINITE F50 PLUS:

Λειτουργία ανακίνησης	Πλάτος ανακίνησης	Συχνότητα ανακίνησης
HIGH (ΥΨΗΛΗ)	2,8 mm	12,3 Hz
NORMAL (KANONIKH)	4,4 mm	9,2 Hz
LOW (XAMH/H)	4,4 mm	7,8 Hz
WIDE (EYPEIA)	14,1 mm	2,0 Hz

4.2 Περιγραφή οργάνου

Η παρακάτω εικόνα παρουσιάζει τα τμήματα του οργάνου.



4. Λειτουργίες οργάνου

τροφοδοσίας



Η φωτεινή ένδειξη (LED) κατάστασης παρέχει πληροφορίες για την κατάσταση του οργάνου:

- Αναβοσβήνει πράσινο: το όργανο δεν είναι συνδεδεμένο με το Magellan •
- Πράσινο: το όργανο είναι συνδεδεμένο και έτοιμο για μέτρηση
- Κόκκινο: μέτρηση σε εξέλιξη

Στην αριστερή πλευρά του οργάνου θα βρείτε τη θύρα USB, το διακόπτη κύριας τροφοδοσίας και την υποδοχή κύριας τροφοδοσίας.



Διακόπτης κύριας τροφοδοσίας

Η ετικέτα με τα στοιχεία του οργάνου βρίσκεται στο κάτω μέρος.

Παράδειγμα πινακίδας στοιχείων



Τα περιεχόμενα της πινακίδας στοιχείων (π.χ. όνομα μοντέλου και αριθμός είδους) μπορεί να ποικίλλουν, ανάλογα με το συγκεκριμένο μοντέλο.

Για επισκόπηση των διάφορων οργάνων για τα οποία ισχύουν οι παρούσες οδηγίες χρήσης, δείτε τη Δήλωση Συμμόρφωσης ("Declaration of Conformity") στην τελευταία σελίδα του παρόντος εντύπου.
4.3 Περιγραφή τροχού φίλτρων

Ο τροχός φίλτρων στο βασικό εξοπλισμό του INFINITE F50 PLUS διαθέτει τέσσερα φίλτρα παρεμβολής στενού φάσματος, τα οποία έχουν σταθερό μήκος κύματος (405, 450, 620 και 492 nm). Μπορείτε να εξοπλίσετε τον τροχό φίλτρων με έως και 8 φίλτρα. Για επιπλέον διαθέσιμα φίλτρα, επικοινωνήστε με τον τοπικό αντιπρόσωπο πωλήσεων της Tecan.

Τα φίλτρα του τυπικού τροχού φίλτρων είναι τοποθετημένα ως εξής:

Θέση φίλτρου	Μήκος κύματος φίλτρου
1	405 nm
2	450 nm
3	620 nm
4	492 nm
5 - 8	κενές θέσεις φίλτρων



Όταν επιλέγεται ένα μήκος κύματος για μια μέτρηση, το συγκεκριμένο φίλτρο μεταφέρεται στη δέσμη φωτός με τη μετακίνηση του τροχού στην αντίστοιχη θέση.



Σημείωση Για περισσότερες πληροφορίες σχετικά με τον ορισμό νέου φίλτρου, ανατρέξτε στο κεφάλαιο 7.5.2 Ορισμός φίλτρων.



4.4 Προδιαγραφές οργάνου

Στους παρακάτω πίνακες αναγράφονται οι προδιαγραφές του μετρητή απορρόφησης INFINITE F50 PLUS.

4.4.1 Γενικές προδιαγραφές

ΠΑΡΑΜΕΤΡΟΙ	ΧΑΡΑΚΤΗΡΙΣΤΙΚΑ
Κύρια τροφοδοσία Εξωτερικό τροφοδοτικό	Παροχή τροφοδοσίας: Βασικό όργανο με προσαρμογέα AC: 100-240 V AC, 50/60 Hz, μέγ. 1,2 A (αυτόματη ανίχνευση τάσης, Κατηγορία υπέρτασης II) Βασικό όργανο χωρίς προσαρμογέα AC: 24 V DC (Κατηγορία υπέρτασης I)
Κατανάλωση INFINITE F50 PLUS	Κατάσταση αναμονής: Περίπου 12 VA Κατάσταση λειτουργίας: Μέγ. 30 VA
Εξωτερικές διαστάσεις	Πλάτος: 34,7 cm(13,66 ίντσες) Βάθος: 18,9 cm(7,44 ίντσες) Ύψος: 13,4 cm(5,28 ίντσες)
Βάρος	2,6 κιλά (συμπεριλαμβάνεται και το τροφοδοτικό)
Θερμοκρασία περιβάλλοντος:	
Λειτουργία	15 °C έως 35 °C (59 °F έως 95 °F)
Αποθήκευση	-30 °C έως 60 °C (-22 °F έως 140 °F)
Σχετική υγρασία:	20% έως 80%
Βαθμός ρύπανσης	2
Μέθοδος απόρριψης	Μολυσμένα απόβλητα
Περιβάλλον	Για περισσότερες πληροφορίες, ανατρέξτε στο κεφάλαιο 3.3 Περιβαλλοντικές απαιτήσεις.



4.4.2 Προδιαγραφές μέτρησης

ΠΑΡΑΜΕΤΡΟΙ	ΧΑΡΑΚΤΗΡΙΣΤΙΚΑ
Χρόνοι μέτρησης: μονό μήκος κύματος διπλό μήκος κύματος	< 15 δευτερόλεπτα < 20 δευτερόλεπτα
Εύρος μήκους κύματος: Τυπικό	400 - 750 nm
Εύρος μέτρησης: 400 - 750 nm	0 – 4,000 OD
Ανάλυση:	0,0001 OD
Ορθότητα: 450, 492 nm 0,000 – 2,000 OD 2,000 – 3,000 OD	≤ (1,0 % + 0,010 OD)* ≤ (1,5 % + 0,010 OD)*
Ακρίβεια: 450, 492 nm 0,000 – 2,000 OD 2,000 – 3,000 OD	≤ (0,5 % + 0,005 OD)* ≤ (1,0 % + 0,005 OD)*
Γραμμικότητα: 450, 492 nm 0,000 – 2,000 OD 2,000 – 3,000 OD	≤ 1 % ≤ 1,5 %
Επιλογή μήκους κύματος: Τυπικό φίλτρο	Φίλτρα παρεμβολής στενού εύρους φάσματος. Μπορούν να τοποθετηθούν έως 8 φίλτρα σε έναν τροχό φίλτρων.
Ορθότητα μήκους κύματος φίλτρων:	Κεντρικό μήκος κύματος ± 2 nm
Εύρος φάσματος φίλτρων: Σε διαπερατότητα 50%	10 ± 2 nm
Πηγή φωτός:	LED
Διασύνδεση υπολογιστή:	USB

Όλες οι συνδεδεμένες συσκευές πρέπει να έχουν εγκριθεί και να καταγράφονται στο IEC 60950-1 Information Technology Equipment – Safety και σε ισοδύναμα τοπικά πρότυπα.

* μεγαλύτερη από ή ίση με το x% της τιμής μέτρησης συν την αντίστοιχη τιμή OD.

4.4.3 Μικροπλάκες

Όλες οι μικροπλάκες 96 κυψελίδων με διάφανο πυθμένα (επίπεδο, σχήματος C, U και V, συμπεριλ. μικροπλακών λωρίδων κυψελίδων) που συμμορφώνονται με τα παρακάτω πρότυπα μπορούν να χρησιμοποιηθούν με το μετρητή απορρόφησης INFINITE F50 PLUS:

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



ΠΡΟΣΟΧΗ

ΧΡΗΣΙΜΟΠΟΙΕΙΤΕ ΜΟΝΟ ΜΙΚΡΟΠΛΑΚΕΣ ΧΩΡΙΣ ΚΑΛΥΜΜΑΤΑ. ΜΗ ΧΡΗΣΙΜΟΠΟΙΕΙΤΕ ΜΙΚΡΟΠΛΑΚΕΣ ΥΨΗΛΟΤΕΡΕΣ ΑΠΟ 15,2 MM.

Χειρισμός της μικροπλάκας

Τοποθετείτε ή αφαιρείτε τη μικροπλάκα μόνο όταν ο φορέας μεταφοράς της πλάκας έχει πλήρως εξαχθεί (όπως εικονίζεται παρακάτω) και ο μηχανισμός μεταφοράς δεν είναι ενεργός. Μην ανοίγετε το κάλυμμα του περιβλήματος ενώ η φωτεινή ένδειξη (LED) κατάστασης είναι κόκκινη.



ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΚΑΤΑ ΤΟ ΧΕΙΡΙΣΜΟ ΤΗΣ ΜΙΚΡΟΠΛΑΚΑΣ, ΝΑ ΦΟΡΑΤΕ ΠΑΝΤΑ ΓΑΝΤΙΑ ΜΙΑΣ ΧΡΗΣΗΣ ΚΑΙ ΠΡΟΣΤΑΤΕΥΤΙΚΟ ΡΟΥΧΙΣΜΟ.



Θέση φρεατίου Α1 Φορέας μεταφοράς πλάκας – σε πλήρη εξαγωγή

4.5 Παρελκόμενα οργάνου

Η παρακάτω λίστα περιλαμβάνει τα διαθέσιμα προαιρετικά παρελκόμενα για το INFINITE F50 PLUS, τα οποία μπορείτε να παραγγείλετε ξεχωριστά:

- Επιπλέον φίλτρα
- Ειδικό εργαλείο τοποθέτησης φίλτρων
- Πλάκα MultiCheck[™] για τη σειρά INFINITE F50

Για περισσότερες πληροφορίες και διαθεσιμότητα στη χώρα σας, επικοινωνήστε με τον τοπικό αντιπρόσωπο πωλήσεων της Tecan.





5. Έλεγχος ποιότητας

5.1 Εισαγωγή



ΠΡΟΣΟΧΗ ΕΑΝ ΟΠΟΙΑΔΗΠΟΤΕ ΣΤΙΓΜΗ, ΥΠΑΡΧΟΥΝ ΑΜΦΙΒΟΛΙΕΣ ΓΙΑ ΤΗΝ ΑΠΟΔΟΣΗ ΑΝΑΛΥΣΗΣ ΤΟΥ INFINITE F50 PLUS, ΑΚΟΛΟΥΘΗΣΤΕ ΤΙΣ ΟΔΗΓΙΕΣ ΠΟΥ ΠΑΡΕΧΟΝΤΑΙ ΓΙΑ ΤΟΝ ΕΛΕΓΧΟ ΠΟΙΟΤΗΤΑΣ Ή ΕΠΙΚΟΙΝΩΝΗΣΤΕ ΜΕ ΤΟ ΤΟΠΙΚΟ ΤΜΗΜΑ ΥΠΟΣΤΗΡΙΞΗΣ ΠΕΛΑΤΩΝ ΤΗΣ ΤΕCAN.

Αυτό το κεφάλαιο παρέχει πληροφορίες σχετικά με τη διαδικασία αυτόματου ελέγχου του οργάνου και οδηγίες για εύκολο έλεγχο της ποιότητας λειτουργίας.

5.2 Διαδικασία αυτόματου ελέγχου

Κατά τη σύνδεση του INFINITE F50 PLUS με το λογισμικό ελέγχου Magellan του μετρητή, ελέγχονται οι μηχανισμοί και οι αισθητήρες και αρχικοποιούνται ο φορέας της πλάκας και ο τροχός φίλτρων.

Πριν από κάθε μέτρηση, εκτελείται μια διαδικασία βαθμονόμησης αυτόματου ελέγχου ώστε να εξασφαλιστεί η σωστή λειτουργία του οργάνου και να βαθμονομηθεί το σύστημα οπτικών στοιχείων.

5.3 Επαλήθευση λειτουργικότητας

Είναι δυνατή η εκτέλεση των παρακάτω ελέγχων για την εξασφάλιση της σωστής λειτουργίας του οργάνου και της λήψης ορθών αποτελεσμάτων.

Η δυνατότητα αναπαραγωγής και ορθότητας των αποτελεσμάτων του οργάνου ενδέχεται να διαφέρει ανάλογα με τον τύπο του διαλύματος και της μικροπλάκας που χρησιμοποιούνται.

Για την εξάλειψη αυτής της επίδρασης, τα όργανα ελέγχονται στο εργοστάσιο με μια πλάκα βαθμονόμησης, η οποία καταργεί την επίδραση του διαλύματος και οποιασδήποτε απόκλισης εξαιτίας της τοποθέτησης της μικροπλάκας κατά τη μέτρηση.

5.3.1 Έλεγχος πολλαπλών σημείων (MultiCheck)

Ο έλεγχος MultiCheck παρέχει έναν αυτοματοποιημένο έλεγχο της απόδοσης του μετρητή, συμπεριλαμβανομένων της ορθότητας, της γραμμικότητας, της ακρίβειας και ευθυγράμμισης με πρότυπα ανιχνεύσιμα στο NIST.

5.3.2 Έλεγχος μικροπλάκας

Εάν οι οπτικές πυκνότητες των κυψελίδων στη μικροπλάκα δεν είναι σταθερές, τα αποτελέσματα που θα ληφθούν με αυτόν τον τύπο μικροπλάκας θα έχουν επηρεαστεί.

Αυτή η έλλειψη σταθερότητας μπορεί να ελεγχθεί με την ανάγνωση μιας κενής μικροπλάκας.

Οι τιμές OD που λαμβάνονται από την κενή μικροπλάκα θα πρέπει να βρίσκονται εντός περιορισμένου εύρους. Για παράδειγμα: ± 0,010 OD.

Εάν οι τιμές OD δεν είναι εντός αυτού του εύρους, αυτός ο τύπος μικροπλάκας δεν πρέπει να χρησιμοποιηθεί.



Με τη χρήση μετρήσεων διπλού μήκους κύματος, η επίδραση της διαφοράς σε τιμές OD της μικροπλάκας καταργείται ή μειώνεται σε βαθμό εντός αποδεκτών ορίων.

5.3.3 Ακρίβεια οργάνου με υγρά δείγματα

Αυτή η διαδικασία μπορεί να χρησιμοποιηθεί για τον έλεγχο της δυνατότητας αναπαραγωγής των μετρήσεων. Συνιστάται η χρήση μικροπλάκας με επίπεδο πυθμένα.

Γεμίστε μια καινούρια μικροπλάκα με ένα μόλις παρασκευασμένο διάλυμα Orange G. Χρησιμοποιήστε διαφορετικές αραιώσεις του διαλύματος σε κάθε κυψελίδα έτσι ώστε να ληφθεί ένα εύρος από οπτικές πυκνότητες. Βεβαιωθείτε ότι οι κυψελίδες περιέχουν τουλάχιστον 200 μl. Η σειρά αραιώσεων θα πρέπει να είναι εντός του εύρους 0,1 έως 3,0 OD. Για την επίτευξη περίπου 3 OD, συνιστάται η χρήση 125 mg.l⁻¹ Orange G (Sigma, Cat. No. O7252).

Ορίστε έναν κύκλο δοκιμών με τη χρήση του φίλτρου 492 nm και, κατόπιν, μετρήστε τη μικροπλάκα τουλάχιστον τρεις φορές.

Για κάθε κυψελίδα, υπολογίστε:

- τη μέση τιμή OD
- τη σταθερή απόκλιση

Παράδειγμα

Αποτελέσματα 0,000 έως 2,000 OD

Η τυπική απόκλιση κάθε κυψελίδας θα πρέπει να είναι εντός (0,5% + 0,005 OD).

Υπολογισμός μέγιστης επιτρεπτής απόκλισης με τη χρήση 1,000 OD ως μέση τιμή OD:

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

Αποτελέσματα 2,001 έως 3,000 OD

Η τυπική απόκλιση κάθε κυψελίδας θα πρέπει να είναι εντός (1,0 % + 0,005 OD).

Υπολογισμός μέγιστης επιτρεπτής απόκλισης με τη χρήση 2,400 OD ως μέση τιμή OD:

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

Αποτελέσματα πάνω από 3,000 OD

Αποτελέσματα πάνω από 3,000 OD χρησιμοποιούνται μόνο ως ένδειξη και η ακρίβεια δεν είναι εγγυημένη.



5.3.4 Γραμμικότητα οργάνου με υγρά δείγματα

Η γραμμικότητα για το όργανο και την εφαρμογή στο μήκος κύματος που χρησιμοποιείται μπορεί να ελεγχθεί με τη χρήση μιας σειράς αραιώσεων ενός διαλύματος.

Το αποτέλεσμα εξαρτάται από την καθαρότητα της χρωστικής που χρησιμοποιείται και του μηνίσκου του υγρού στις κυψελίδες.

Ως αναφορά μπορεί να χρησιμοποιηθεί μια σειρά αραιώσεων του διαλύματος Orange G για μετρήσεις στα 492 nm.

Η σειρά αραιώσεων θα πρέπει να είναι εντός του εύρους 0,1 έως 3,0 OD. Για την επίτευξη περίπου 3 OD, συνιστάται η χρήση 125 mg.l⁻¹ Orange G (Sigma, Cat. No. 07252).

Για άλλα μήκη κύματος, πρέπει να χρησιμοποιηθούν διαφορετικά διαλύματα.

Κατόπιν, 200 μΙ από κάθε αραίωση μεταφέρονται με πιπέτα στη μικροπλάκα. Θα πρέπει να χρησιμοποιηθούν τουλάχιστον δύο δείγματα από κάθε αραίωση για τη μείωση των σφαλμάτων που προκαλούνται από τη χρήση πιπέτας.

Ακολουθεί η μέτρηση της μικροπλάκας και εξάγεται η γραμμική παλινδρόμηση OD σε σχέση με τη συγκέντρωση, από το μέσο όρο των μετρημένων τιμών OD.

Καθορίστε τη μη σταθμισμένη τιμή τετραγώνου υπολοίπου R² της γραμμής παλινδρόμησης.

Οι τυπικές τιμές τετραγώνου υπολοίπου για μια τυπική εφαρμογή είναι ίσες με ή μεγαλύτερες από R² = 0,998.



Σημείωση Τα δεδομένα μπορεί να διαφέρουν εξαιτίας ανακριβειών από τη χρήση πιπέτας.



6.1 Εισαγωγή

Τα **αρχεία παραδειγμάτων** του Magellan προσφέρουν μεθόδους και χώρους εργασίας Magellan κάνοντας μια εισαγωγή στο λογισμικό και διευκολύνοντας το χρήστη στην εργασία με αυτό. Τα αρχεία παραδειγμάτων για έναν ποσοτικό και ποιοτικό προσδιορισμό ELISA εγκαθίστανται αυτόματα με την εγκατάσταση του Magellan.

6.2 Παράδειγμα «βήμα-προς-βήμα»: Ποσοτικός προσδιορισμός ELISA

Σε αυτό το κεφάλαιο παρέχεται ένα παράδειγμα «βήμα-προς-βήμα» (ποσοτικός έλεγχος) για τον τρόπο δημιουργίας μεθόδου στο Magellan. Ακολουθώντας τις οδηγίες θα μάθετε πώς να ορίζετε αξιολογήσεις από περιγραφή του κιτ ελέγχου στο Magellan.



Σημείωση Τα αρχεία παραδειγμάτων εμφανίζονται αυτόματα στο Method List (Λίστα μεθόδων) στο Magellan. Στο Magellan Tracker, αυτά τα αρχεία είναι διαθέσιμα στην προεπιλεγμένη διαδρομή δεδομένων και πρέπει να μετατραπούν.

6.2.1 Περιγραφή κιτ ελέγχου

Στην περιγραφή του κιτ ελέγχου του κατασκευαστή για ένα ποσοτικό προσδιορισμό ELISA ανίχνευσης αντισωμάτων IgM, υπάρχουν οι παρακάτω οδηγίες: Διάταξη πλάκας

	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										
		1	1	1	1	1	1	1	1	1	1	t

BLK = Τυφλές, NC = Αρνητικός μάρτυρας, C1 - C6 = Βαθμονομητές (Πρότυπα), S1 - S… = Δείγματα



Μέτρηση και αξιολόγηση

Πραγματοποιήστε μέτρηση της πλάκας σε μήκος κύματος 492 nm, με αναφορά στα 620 nm.

Μηδενίστε το μετρητή/την πλάκα στην κυψελίδα Α1.

Συγκεντρώσεις βαθμονομητών (Πρότυπα):

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

Μετά τη διόρθωση τυφλού, οι οπτικές πυκνότητες (OD 492 - OD 620) σχεδιογραφούνται σε σχέση με τη συγκέντρωση. Η γραμμή παλινδρόμησης που διαπερνά αυτά τα σημεία είναι η πρότυπη καμπύλη.

Ερμηνεία των αποτελεσμάτων ελέγχου:

lgM < 18 UA/ml	Αρνητικό
18 UA/ml ≤ IgM < 22 UA/ml	Ενδιάμεσο
IgM ≥ 22 UA/mI	Θετικό

Η υπολογισμένη συγκέντρωση IgM και των δύο αρνητικών μαρτύρων πρέπει να είναι κάτω από 8 UA/ml.

Χειρισμός δεδομένων

Μετά τη μέτρηση, το αρχείο δεδομένων (χώρος εργασίας) αποθηκεύεται αυτόματα και δημιουργείται μια αναφορά που περιέχει τις παραμέτρους μέτρησης, τη διάταξη πλάκας, τις τιμές τυφλών, την πρότυπη καμπύλη, τις συγκεντρώσεις IgM, τον ορισμό ορίων, τα ποιοτικά αποτελέσματα των δειγμάτων και των επικυρώσεων.

Επιπλέον, η διάταξη και τα ποιοτικά αποτελέσματα αποθηκεύονται ως αρχείο ASCII.



6.2.2 Δημιουργία μεθόδου

Στο πλαίσιο διαλόγου Wizard List (Λίστα οδηγών), επιλέξτε Create/edit a method (Δημιουργία/επεξεργασία μεθόδου) και κάντε κλικ στο OK. Κάντε κλικ στο Continue (Συνέχεια) στη σελίδα Welcome (Καλώς ορίσατε) στον οδηγό Create/edit a method (Δημιουργία/επεξεργασία μεθόδου). Εμφανίζεται το πλαίσιο διαλόγου Select a file (Επιλογή αρχείου). Επιλέξτε New (Δημιουργία).

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

Παράμετροι μέτρησης

Κάντε κλικ στο **Make Your Selection** (Επιλογή). Εμφανίζεται η σελίδα **Measurement parameters** (Παράμετροι μέτρησης).





Στη λωρίδα **Wavelength** (Μήκος κύματος), επιλέξτε 492 nm ως μήκος κύματος στο πεδίο Measurement (Μέτρηση) και 620 nm ως μήκος κύματος στο πεδίο Reference (Αναφορά).

🖌 🔻 Absorb	ance	3	}
-Wavelength-		Label	
Measurement:	492 nm 🖌 🖌	Name: Label1	
Reference:	620 nm 🔽		

Συνεχίστε τον Οδηγό κάνοντας κλικ στο **Choose measurement parameters** (Επιλογή παραμέτρων μέτρησης). Εμφανίζεται το παράθυρο **Plate layout** (Διάταξη πλάκας).





Σχεδίαση διάταξης

Ορίστε τη διάταξη της πλάκας με τη χρήση του πλαισίου διαλόγου **Well Assignment** (Αντιστοίχιση κυψελίδων) στη δεξιά πλευρά της οθόνης.

We	ell Assig	ınm.	
C	Identifier	s	Exp. group:
	SM	^	1
	BL		-ID-Num.
	ST		≠ 4
	PC		1
	NC		
	LPC		- Replicates -
	HPC		≠ 4
	BF		Ix number:
	RF	~	1
I	Def. ide	ntif.	O All
	Fill selecti	on	Delete selection

Στο πλαίσιο ομάδας Identifiers (Αναγνωριστικά), επιλέξτε BL (Blank) (Τυφλό).

Στο πλαίσιο ομάδας Experimental (Πειραματικές) παραμένει ο αριθμός 1.

Στο πλαίσιο ομάδας **Replicates** (Αντίγραφα), η επιλογή **All** (Όλες) έχει οριστεί αυτόματα.

Κάντε κλικ στην κυψελίδα **Α1**, η οποία στη συνέχεια επισημαίνεται με ένα κόκκινο πλαίσιο.

Κάντε κλικ στο **Fill selection** (Πλήρωση με επιλεγμένο) και η κυψελίδα παίρνει το όνομα του επιλεγμένου τύπου αναγνωριστικού.



Σημείωση Μια κυψελίδα μπορεί επίσης να πληρωθεί με διπλό κλικ σε αυτή.

Επιλέξτε τις παρακάτω ρυθμίσεις στο πλαίσιο διαλόγου Well Assignment (Αντιστοίχιση κυψελίδων):

Στο πλαίσιο ομάδας **Identifiers** (Αναγνωριστικά), επιλέξτε **NC** (Negative Control) (Αρνητικός μάρτυρας).

Στο πλαίσιο ομάδας Experimental (Πειραματικές) παραμένει ο αριθμός 1.

Στο πλαίσιο ομάδας **Replicates** (Αντίγραφα), η επιλογή **All** (Όλες) έχει οριστεί αυτόματα.

Ξεκινώντας από την κυψελίδα **B1** κάντε κλικ και σύρετε το ποντίκι στη **C1**. Οι κυψελίδες **B1** έως **C1** επισημαίνονται με ένα κόκκινο πλαίσιο.

Κάντε κλικ στο **Fill selection** (Πλήρωση με επιλεγμένο) και οι κυψελίδες επισημαίνονται με τον επιλεγμένο τύπο αναγνωριστικού.

Στις κυψελίδες **D1** έως **G2** πρέπει να αντιστοιχιστούν βαθμονομητές (τυπικοί). Επιλέξτε τις παρακάτω ρυθμίσεις στο πλαίσιο διαλόγου **Well Assignment** (Αντιστοίχιση κυψελίδων):



Στο πλαίσιο ομάδας **Identifiers** (Αναγνωριστικά), επιλέξτε **ST (Standard)** (Πρότυπα).

Στο πλαίσιο ομάδας Experimental (Πειραματικές) παραμένει ο αριθμός 1.

Στο πλαίσιο ομάδας **Replicates** (Αντίγραφα), επιλέξτε μεταξύ **Fix number** (Συγκεκριμένος αριθμός) και **All** (Όλες).

Fix number (Συγκεκριμένος αριθμός):

Ενεργό μόνο για πρότυπα και δείγματα όπου μπορούν να χρησιμοποιηθούν αναγνωριστικά.

Εάν το κουμπί **Fix number** (Συγκεκριμένος αριθμός) είναι ενεργό, μπορείτε να εισαγάγετε έναν αριθμό στο αντίστοιχο πεδίο κειμένου. Αυτός ο αριθμός καθορίζει πόσα αντίγραφα προορίζονται για αυτήν τη μέθοδο. Στις επιλεγμένες κυψελίδες, δημιουργείται για κάθε αναγνωριστικό ο αριθμός αντιγράφων που εισήγατε. Κατά συνέπεια, ο αριθμός των επιλεγμένων κυψελίδων πρέπει να είναι πολλαπλάσιο του αριθμού αντιγράφων που εισήγατε.

All (Όλες):

Όλες οι επιλεγμένες κυψελίδες ορίζονται ως αντίγραφα. Εάν έχει επιλεγεί ένας υπάρχων αριθμός αναγνωριστικού για τα δείγματα και τα πρότυπα, οι επιλεγμένες κυψελίδες προστίθενται ως αντίγραφα στα υπάρχοντα αντίγραφα. Με όλους τους άλλους τύπους αναγνωριστικών, οι επιλεγμένες κυψελίδες προστίθενται ως αντίγραφα στα υπάρχοντα αντίγραφα.

Δύο κουμπιά με βελάκια σρίζουν την ακολουθία των αντιγράφων και των αριθμών αναγνωριστικών (οριζόντια ή κάθετα).

Σε αυτό το παράδειγμα, επιλέξτε Fix Number (Συγκεκριμένος αριθμός) και 2.

Στο πλαίσιο **ID-Number** (Αριθμός αναγνωριστικών) και στο πλαίσιο ομάδας **Replicates** (Αντίγραφα), επιλέξτε τα **κάθετα βέλη**.

Κατόπιν, επιλέξτε τις κυψελίδες D1 έως G2 και κάντε κλικ στο **Fill selection** (Πλήρωση με επιλεγμένο).



Σημείωση

Επιλέξτε τις κυψελίδες ως εξής: Ξεκινώντας από την κυψελίδα D1, κάντε κλικ και σύρετε το ποντίκι πάνω από τις απαιτούμενες κυψελίδες έως την H1. Κατόπιν, κρατήστε πατημένο το πλήκτρο Ctrl και σύρετε το ποντίκι πάνω από τις απαιτούμενες κυψελίδες από την A2 έως την G2.



Create/Edit a Method															
Undo Redo Select al unu	sed	10%	100% Zo	om											
(Ś														
Method layout 🔗	5														
Plate layout	ate/														
Conc, Dil, Refvalues	edit		-	0	0		P		-	0		10	44	40	$\setminus \parallel$
粒 Transformed data 🛞	meth		1	2	3	4	5	0		8	g	10	11	12	
Add new transformation	8		DI 4	074 0									Well Assign	n.	
Concentrations 🔗	ΓII		1/1	2/2									Identifiers	Exp. group:	
Standard curve		0-0											SM S		
Evaluate data 🔗			NC1	ST1_4									ST	ID-Num. ⇒+ Ia	
Cutoff definition			1/2										PC	7	
			NOI	074 4									NC		
Data handling 🔗		C	NC1 2/2	2/2									LPC	Replicates	
Printed report			0.2										RF	44	
Automated data handling			ST1_1	ST1_5									RF	2	я.
🧭 Miscellaneous 🔗			1/2										Def identi		
User prompts Number format															
Method notes		E	ST1_1	ST1_5									Fill	Delete	
		5	212												-'
			ST1 2	ST1 6									i		
		F	1/2	1/2											
		- T													
			ST1_2	ST1_6											Istria
		G	2/2	2/2											N AL
			CT1 2											_	CAL
		H	1/2												μ
		00													
															/ J
		_			_			_					_		
Help														NEXT	
Cancel <<< Back															
													@ 2005	Tecan	

Εμφανίζεται η σελίδα Plate Layout (Διάταξη πλάκας), ως εξής:

Κάντε κλικ στο **Select all unused** (Επιλογή όλων των μη χρησιμοποιούμενων) από τη γραμμή εργαλείων για να επιλέξετε όλες τις κενές κυψελίδες στην πλάκα. Κατόπιν, κρατήστε πατημένο το πλήκτρο Ctrl και κάντε κλικ στην κυψελίδα **H12**, έτσι ώστε να παραμείνει κενή και χωρίς επισήμανση.

Στο πλαίσιο διαλόγου **Well Assignment** (Αντιστοίχιση κυψελίδων), επιλέξτε **SM** (Sample) (Δείγμα) κάτω από το *Identifiers (Αναγνωριστικά*).

Στο πλαίσιο ομάδας Experimental (Πειραματικές) παραμένει ο αριθμός 1.

Στο πλαίσιο ομάδας **Replicates** (Αντίγραφα), επιλέξτε μεταξύ **Fix number** (Συγκεκριμένος αριθμός) και **2**.

Στο πλαίσιο **ID-Number** (Αριθμός αναγνωριστικών) αφήστε το 1 και στο πλαίσιο ομάδας **Replicates** (Αντίγραφα), επιλέξτε τα **κάθετα βέλη**. Κάντε κλικ στο **Fill selection** (Πλήρωση με επιλεγμένο). Η διαδικασία ορισμού διάταξης έχει ολοκληρωθεί.



Μετασχηματισμοί

Στη γραμμή ελέγχου, στα αριστερά του παραθύρου, προχωρήστε στην επόμενη επιλογή, Add new transformation... (Προσθήκη νέου μετασχηματισμού) από το στοιχείο Transformed data (Μετασχηματισμένα δεδομένα), ώστε να ορίσετε μια αναγωγή της απορρόφησης βάσει τυφλών.

Εμφανίζεται πλαίσιο διαλόγου που σας ρωτά εάν θέλετε να ορίσετε μια αναγωγή της απορρόφησης βάσει των τυφλών. Κάντε κλικ στο **Yes** (Nai). Εμφανίζεται το παρακάτω παράθυρο:

Create/Edit a Method															
Undo Redo		10%	100% Zo	om											
S		Input data:	Difference data				~							Constants	Options
Method layout 🛞	8	fx	×8L1							~	Available dat	a (multiple data	sets) 🔻	Functions&Co	nstants 🔻
Plate layout	ate/e														
Conc.r, bitr, Keitrvaldes	dit n														
Blank reduction	etho		പ	9	2	4	E	R	7	0	0	40	ଶାଶ	สด	
Add new transformation	7		U	6	3	43	2	0	Ű	0	8		0.0	12)
Concentrations			BL1	ST1 3	SM1 1	SM1 5	SM1 9	SM1 13	SM1 17	SM1 21	SM1 25	SM1 29	SM1 33	SM1 37	
Standard curve		A	1/1	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Evaluate data 🔗			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Cutoff definition			NC1	ST1_4	SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
QC Valdation			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Data handling 🛞			NC1	ST1_4	SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
Printed report			2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Automated data handling			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
🧱 Miscellaneous 🛞			SI1_1 1/2	SI1_5 1/2	SM1_3	SM1_7 1/2	SM1_11 1/2	SM1_15 1/2	SM1_19 1/2	SM1_23	SM1_27 1/2	SM1_31	SM1_35	SM1_39 1/2	
User prompts Number format			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Method notes			ST1_1	ST1_5	SM1_3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
			2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
			X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	X-BL1	
		F	1/2	511_6 1/2	5MI1_4 1/2	5MI_8 1/2	SM1_12 1/2	SM1_10 1/2	5MI1_20	SM1_24	SM1_28	5M1_32 1/2	5M1_30 1/2	5M1_40 1/2	
		0	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
			ST1_2	ST1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	stria
		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	4 Au
			X-BLI	X-BLT	X-BLI	X-BLT	X-BLI	X-BLI	X-BLT	X-BLI	X-BLI	X-BLI	X-BLI	X-BL1	CAN
		H H	1/2	5MIT_1 1/2	5IVIT_5	5MT_9	1/2	5WI1_17 1/2	1/2	5MT_25 1/2	5MT_29 1/2	1/2	1/2		μ
		00	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1		
Help															
Cancel (<< Back														NEXT	

Το Difference Data (Δεδομένα διαφοράς) έχει επιλεγεί αυτόματα στο πλαίσιο **Input data** (Δεδομένα εισόδου). Εάν έχετε επιβεβαιώσει τον ορισμό μιας μείωσης τυφλών προηγουμένως, το λογισμικό αυτόματα το ονομάζει **Blank reduction** (Αναγωγή απορρόφησης βάσει τυφλών) (δείτε τα μετασχηματισμένα δεδομένα στη γραμμή ελέγχου).

Στο πλαίσιο **Formula** (Τύπος) εμφανίζεται αυτόματα το **x-BL1** για αυτή την αναγωγή απορρόφησης βάσει των τυφλών, όπου x είναι η τρέχουσα τιμή δεδομένων εισόδου σε μια κυψελίδα και BL1 είναι η μέση τιμή των τυφλών κυψελίδων της πειραματικής ομάδας 1.

Για περισσότερες λεπτομέρειες και επεξηγήσεις σχετικά με τον ορισμό και την αντιστοίχιση μετασχηματισμών, ανατρέξτε τις Οδηγίες Χρήσης του Magellan.

Σε κάθε κυψελίδα εμφανίζονται οι παρακάτω πληροφορίες (παράδειγμα κυψελίδας A5):

SM1_9	Δείγμα, αριθμός πειραματικής ομάδας 1, αριθμός αναγνωριστικού δείγματος 9.
2/2	Αριθμός αντιγράφου 2, συνολικός αριθμός αντιγράφων 2.
x-BL1 ή 1	Αντιστοιχισμένος μετασχηματισμός x-BL1 (όταν έχει επιλεγεί Transformation (Μετασχηματισμός) ή τιμή Dilution Factor (Παράγων αραίωσης) 1 (όταν έχουν επιλεγεί τιμές Conc., Dil,. Ref. (Συγκέντρωση, Αραίωση, Αναφορά)).



Ορισμός τιμών Συγκέντρωσης / Αραίωσης / Αναφοράς

Στη γραμμή ελέγχου, επιλέξτε **Conc., Dil., Ref.-values** (Τιμές Συγκέντρωσης, Αραίωσης, Αναφοράς) από το στοιχείο **Method layout** (Διάταξη μεθόδου) για να ορίσετε τις αντίστοιχες τιμές, όπως περιγράφεται στο κιτ ελέγχου.

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

Βεβαιωθείτε ότι έχει επιλεγεί το **ST** (Πρότυπα) στη λίστα **Select Identifier** (Επιλογή αναγνωριστικού).

Στη λίστα **Identifier** (Αναγνωριστικά), εμφανίζεται μια λίστα προτύπων από την πειραματική ομάδα 1. Στο αντίστοιχο πλαίσιο **Concentration** (Συγκέντρωση) του **ST1_1** πληκτρολογήστε τον αριθμό **5** και στο πλαίσιο **Unit** (Movάδα) πληκτρολογήστε UA/ml. Στο αντίστοιχο πλαίσιο **Concentration** (Συγκέντρωση) του **ST1_2** πληκτρολογήστε τον αριθμό **10**. Ο ορισμός της μονάδας πρέπει να γίνει μόνο μία φορά και ισχύει για όλα τα πρότυπα. Παρομοίως, πληκτρολογήστε τις τιμές για τα ST1_3 έως ST1_6.

Εμφανίζεται η οθόνη που απεικονίζει τη διάταξη της πλάκας και τη συγκέντρωση.



Πρότυπη καμπύλη

Στη γραμμή ελέγχου, κάντε κλικ στο **Standard curve** (Πρότυπη καμπύλη) από το στοιχείο **Concentrations** (Συγκεντρώσεις) για να ορίσετε την κατάλληλη πρότυπη καμπύλη.

Τα παρακάτω βρίσκονται στην περιγραφή του κιτ ελέγχου:

Μετά τη διόρθωση βάσει τυφλών, οι οπτικές πυκνότητες (OD 492 - OD 620) σχεδιογραφούνται σε σχέση με τη συγκέντρωση. Η γραμμή παλινδρόμησης που διαπερνά αυτά τα σημεία είναι η πρότυπη καμπύλη.



Στην καρτέλα **Data** (Δεδομένα), επιλέξτε **Blank reduction** (Αναγωγή απορρόφησης βάσει τυφλών) ως δεδομένα εισόδου.

Data	Analysis type Intercepts Axis Graph
	Input data: Blank reduction
	Standards from ext. file: Select
	Standards from experimental group:
	No Standard curve graph
	Additional concentrations

Στην καρτέλα **Analysis type** (Τύπος ανάλυσης), επιλέξτε Linear regression (Γραμμική παλινδρόμηση).

Analysis Type Point to point Linear regression Non-linear regression Cubic spline Akima	Data scaling: Lin(x)Lin(y) 🗸
 Polygomial Four parameters Four parameters Marquardt Five parameters LogitLog 	More
Include (0,0) Extrapolation factor:	



Στην καρτέλα **Axis** (Άξονας), ορίστε τη σήμανση και την κλιμάκωση του άξονα, όπως εικονίζεται παρακάτω:

Label: Concentration [UA/m] Color:	
Color: Log. Scaling O Auto select range Min.: Max.: Max.:	
Auto select range Range Min.: Max.:	
O Range Min.: Max.:	
V Grid Color: Style: ✓	
^-axis	
Label: Blank reduction	
Color: Log. Scaling	
• Auto select range	
O Range Min.: Max.:	

Στην καρτέλα **Graph** (Γράφημα), ορίστε τον τίτλο του γραφήματος, τις καμπύλες, τη γραμματοσειρά και τι θα εμφανίζεται με το γράφημα.

Color:			
Curves			
	Label: Grp. 1		
Color:	Hide curve		
Symbol:	Line width: 1		
Size: Medium 💌			
Font	Display		
⊙ Small	🔽 Legend	✓ Intercepts	
U Medium	🗹 Base points	Error bars	



Ορισμός ορίων

Στη γραμμή ελέγχου, επιλέξτε **Cutoff definition** (Ορισμός ορίων) από το στοιχείο **Evaluate data** (Αξιολόγηση δεδομένων) για να ορίσετε τα όρια της ποιοτικής αξιολόγησης.

Η περιγραφή του κιτ ελέγχου περιλαμβάνει τις παρακάτω οδηγίες:

Ερμηνεία των αποτελεσμάτων ελέγχου:

IgM < 18 UA/ml	Αρνητικό
18 UA/ml ≤ IgM < 22 UA/ml	Ενδιάμεσο
IgM ≥ 22 UA/mI	Θετικό

Για να ορίσετε τα κατάλληλα όρια, ακολουθήστε την παρακάτω διαδικασία:

Στο πλαίσιο Input data (Δεδομένα εισόδου), επιλέξτε **Mean conc. (UA/mI)** (Μέση συγκέντρωση [UA/mI]).

Ο πίνακας **Cutoffs** (Όρια) αντιπροσωπεύει μια κλίμακα που υποδεικνύει το άνω και το κάτω άκρο για τις επιλογές **Limits** (Όρια) και **Labels** (Σημάνσεις). Στο πεδίο **Limits** (Όρια) πληκτρολογήστε 22 ως το πρώτο (υψηλότερο) όριο και 12 ως το δεύτερο (χαμηλότερο) όριο.

Στο πεδίο Labels (Σημάνσεις), εισαγάγετε την ερμηνεία του ελέγχου (Positive, Intermediate και Negative) (Θετική, Ενδιάμεση και Αρνητική) σε κάθε πλαίσιο. Χρησιμοποιήστε την αναπτυσσόμενη παλέτα χρωμάτων για να αντιστοιχίσετε ένα χρώμα:

- Θετική Κόκκινο
- Ενδιάμεση Μπλε
- Αρνητική Πράσινο

Η οθόνη περιέχει τα παρακάτω:

Colors	Labels	Limits		
-	positive	- 22		
-	intermedial	10		
-	negative			
-			 	
-		-	 	
-		-		
		-		
		-	 	
		-	 	
Formula input				

Κάντε κλικ στο **Cutoff results selection** (Επιλογή αποτελεσμάτων ορίων) για να επιλέξετε τους τύπους αναγνωριστικών για τους οποίους τα αποτελέσματα ορίων πρέπει να εμφανίζονται.



Ορισμός επικυρώσεων ελέγχου ποιότητας

Στη γραμμή ελέγχου, κάντε κλικ στο **QC Validations** (Επικυρώσεις ελέγχου ποιότητας) από το στοιχείο **Evaluate data** (Αξιολόγηση δεδομένων). Τα κριτήρια επικύρωσης για τον έλεγχο θα πρέπει να οριστούν έτσι ώστε η εγκυρότητα των αποτελεσμάτων του ελέγχου να είναι εγγυημένη.

Σε αυτό το παράδειγμα, θα πρέπει να πληρείται η παρακάτω απαίτηση:

Η υπολογισμένη συγκέντρωση IgM και των δύο αρνητικών μαρτύρων πρέπει να είναι κάτω από 8 UA/ml.

Στο πλαίσιο Input data (Δεδομένα εισόδου), επιλέξτε **Single conc. (UA/mI)** (Μία συγκέντρωση [UA/mI]).

Στην πρώτη σειρά, πληκτρολογήστε NC1_1<8



Σημείωση NC1_1 σημαίνει Αρνητικός μάρτυρας πειραματικής ομάδας 1, αντίγραφο 1

Στη δεύτερη σειρά, πληκτρολογήστε NC1_2<8

Το πλαίσιο διαλόγου **QC Validations** (Επικυρώσεις ελέγχου ποιότητας) εμφανίζεται τώρα ως εξής:

alidation Condi	itions			
IC1_1<8				
IC1_2<8				
nula input				
Variable	Operators	E		
	LINCIGUNA	EURCOORS		
BL1	V +	and	~	
BL1	· · ·	and	~	
BL1 e to Plate QC	• +	and	~	
BL1 e to Plate QC		and	~	
BL1 e to Plate QC Input data:	Difference data	and	~	
BL1 to Plate QC Input data: Control 1:	Difference data	And		1
BL1 e to Plate QC Input data: Control 1: Control 2:	Difference data	Mean:		<u>Ť</u>
BL1 a to Plate QC Input data: Control 1: Control 2: Control 2:	Difference data	Mean:		
BL1 a to Plate QC Input data: Control 1: Control 2: Control 3:	Difference data	Mean:		<u>†</u>
BL1 e to Plate QC Input data: Control 1: Control 2: Control 3: Control 4:	Difference data BL1 BL1 BL1 BL1	Mean:		↑
BL1 b to Plate QC Input data: Control 1: Control 2: Control 3: Control 4:	Difference data	Mean:		



6.2.3 Οργάνωση έντυπης αναφοράς

Στη γραμμή ελέγχου, κάντε κλικ στο **Printed repor**t (Έντυπη αναφορά) από το στοιχείο **Data handling** (Χειρισμός δεδομένων). Εμφανίζεται η παρακάτω οθόνη:

O List ○ Matrix	You can use drag & drop to create a report. Pick an item from the available data list and drag it into the selected data list. Drop it into the appropriate position inside the report.			
vailable data:	Selected data:			
	Insert E Data Collection Up Append Image: Collection in the summer parameters Down Image: Collection in the summer parameters Image: Collection in the summer parameters Down Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Down Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Down Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summer parameters Image: Collection in the summe			
E Signature	Import			

Στην καρτέλα **Data selection** (Επιλογή δεδομένων), όλα τα διαθέσιμα δεδομένα αναφορών περιέχονται στο πλαίσιο **Available data** (Διαθέσιμα δεδομένα). Χρησιμοποιώντας τα κουμπιά **Insert** (Εισαγωγή) και **Append** (Επισύναψη), μπορείτε να μεταφέρετε δεδομένα στο πλαίσιο **Selected data** (Επιλεγμένα δεδομένα). Η μεταφορά δεδομένων μπορεί επίσης να γίνει με μεταφορά και απόθεση.

Στο πλαίσιο διαλόγου **Print as** (Εκτύπωση ως), επιλέξτε μεταξύ της εκτύπωσης των δεδομένων ως πίνακα ή ως λίστας με ειδικό προσανατολισμό.

Σε αυτό το παράδειγμα, θα δημιουργηθεί μια αναφορά που θα περιέχει τις παραμέτρους μέτρησης, τη διάταξη πλάκας, τις τιμές τυφλών, την πρότυπη καμπύλη, συγκεντρώσεις IgM, τον ορισμό ορίων, ποιοτικά αποτελέσματα δειγμάτων και επικυρώσεις.

Πριν τη δημιουργία της αναφοράς, η προεπιλογή Vertical list/Difference data (Κάθετη λίστα/Δεδομένα διαφοράς) θα πρέπει να καταργηθεί από το πλαίσιο Selected data (Επιλεγμένα δεδομένα). Μόνο το Measurement parameters (Παράμετροι μέτρησης) παραμένει στον πλαίσιο Selected data (Επιλεγμένα δεδομένα). Η επιλογή Print as List (Εκτύπωση ως λίστα) πρέπει να αλλαχθεί σε Print as Matrix (Εκτύπωση ως πίνακας).



C List Matrix	You can use drag & drop to create a report. Pick an item from the available data list and drag it into the selected data list. Drop it into the appropriate position inside the report.
wailable data:	Selected data:
	Inset Append
E Audit Trail	Import

Επιλέξτε Method layout/Layout (Διάταξη μεθόδου/Διάταξη) στο πλαίσιο Available data (Διαθέσιμα δεδομένα) και επισυνάψτε το ως πίνακα στην αναφορά κάνοντας κλικ στο Append (Επισύναψη) Κατόπιν, εισαγάγετε Blank reduction (Αναγωγή απορρόφησης βάσει τυφλών) Mean conc. (UA/ml) (Μέση συγκέντρωση) και Cutoff results (Αποτελέσματα ορίων) στον πίνακα επιλέγοντας τα αντίστοιχα στοιχεία και κάνοντας κλικ στο Insert (Εισαγωγή).

Χρησιμοποιήστε την επιλογή **Append** (Επισύναψη) για να μεταφέρετε τα **Graph: Standard curve** (Γράφημα: Πρότυπη καμπύλη), **Cutoff definition** (Ορισμός ορίων) και **QC Validation criteria** (Κριτήρια επικύρωσης ελέγχου ποιότητας) στα επιλεγμένα δεδομένα. Το τμήμα ρύθμισης δεδομένων για τη διαδικασία ορισμού της αναφοράς έχει ολοκληρωθεί. Το πλαίσιο διαλόγου **Printed Report** (Έντυπη αναφορά) εμφανίζεται όπως παρακάτω:



Στις καρτέλες **Header** (Κεφαλίδα) και **Footer** (Υποσέλιδο), ορίστε τη διάταξη της κεφαλίδας και του υποσέλιδου της αναφοράς (για περισσότερες πληροφορίες, δείτε τις Οδηγίες Χρήσης Magellan).



Εξαγωγή δεδομένων

Στη γραμμή ελέγχου, επιλέξτε **Data export** (Εξαγωγή δεδομένων) από το στοιχείο **Data handling** (Χειρισμός δεδομένων). Σε αυτό το παράδειγμα, η διάταξη και τα αποτελέσματα ορίων αποθηκεύονται ως αρχείο ASCII. Επιλέξτε **Layout** (Διάταξη) και **Cutoff results** (Αποτελέσματα ορίων) από το παράθυρο **Available data** (Διαθέσιμα δεδομένα). Κάντε κλικ στο βέλος → για να τα εισαγάγετε στο παράθυρο **Selected data** (Επιλεγμένα δεδομένα). Στην οθόνη εμφανίζονται οι παρακάτω πληροφορίες:

Available data:		Selected data:	
🕀 🛍 Instrument data	^	Layout	Up
🖶 🤆 Reduced data		Difference data	Doun
Difference data			DOW
🗐 Difference data - Mean			
📃 Difference data - Standard deviati	10		
Difference data - Variation coeffici	ы		
🗄 🔏 Transformed data			
🖶 🚺 Concentrations		→	
Cutoff results			
∎ ¹ ², Sample IDs		< <u>.</u>	
🖃 🛲 Method layout			
Well positions			
Strip method names			
🔄 Original Concentrations			
Dilution factors	~		
Export Options Export to ASCIL File	Г	Export to Excel	
		Enport to Endot	



Σημείωση Τα δεδομένα που εξάγονται θα πρέπει πάντα να περιέχουν τη διάταξη ή τη λίστα αναγνωριστικών δειγμάτων.



Αυτοματοποιημένος χειρισμός δεδομένων

Στη γραμμή ελέγχου, επιλέξτε Automated data handling (Αυτοματοποιημένος χειρισμός δεδομένων) από το στοιχείο Data handling (Χειρισμός δεδομένων).

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

Επιλέξτε export to ASCII file (εξαγωγή σε αρχείο ASCII) και view results after measurements (προβολή αποτελεσμάτων μετά τις μετρήσεις). Στο Magellan Tracker, η επιλογή save workspace (αποθήκευση χώρου εργασίας) είναι επιλεγμένη από προεπιλογή και δεν μπορεί να τροποποιηθεί.

Αποθήκευση μεθόδου

Κάντε κλικ στο **Next** (Επόμενο) για να ανοίξετε το παράθυρο **Save as** (Αποθήκευση ως). Εισαγάγετε το όνομα αρχείου της μεθόδου και συμπληρώστε οποιοδήποτε άλλο πεδίο χρειάζεται.

Create/Edit a Method							
Save in:	🗁 mth	Name 🔺	Remarks	Status			
Filename:	Method1.mth						
File remark	s:						
Audit trail c	omment:						
Org	anize favorites Audit trail	Signatures	Method password				
H Ca	Help Cancel Cancel Cancel						
			C	2009 Tecan			



Πεδίο κειμένου Filename (Όνομα αρχείου)	Πρέπει να εισαγάγετε ένα όνομα αρχείου. Το πρόγραμμα σας προτείνει αυτόματα ένα προεπιλεγμένο όνομα αρχείου αλλά μπορείτε να το αλλάξετε.
Πεδίο κειμένου File remarks (Σχόλια αρχείου)	Τα σχόλια που εισάγονται εδώ θα αποθηκευτούν και θα εμφανίζονται με το όνομα αρχείου.
Πεδίο κειμένου Audit trail comment (Σχόλιο αναφοράς κινήσεων)	Τα σχόλια που εισάγονται εδώ θα αποθηκεύονται στο αρχείο αναφοράς κινήσεων. Αυτή η λειτουργία είναι διαθέσιμη μόνο στο Magellan Tracker .
Κουμπί Organize Favorites… (Οργάνωση προτιμήσεων…)	Εμφανίζεται το πλαίσιο διαλόγου Organize Favorites (Οργάνωση προτιμήσεων).
Πεδίο κειμένου Method Password (Κωδικός μεθόδου)	Εισαγάγετε έναν κωδικό μεθόδου για να προστατέψετε τη μέθοδο.
Πλαίσιο ελέγχου Run this method now (Εκτέλεση αυτής της μεθόδου τώρα)	Αυτή η μέθοδος θα εκτελεστεί αμέσως αφού κάνετε κλικ στο Save&Finish (Αποθήκευση και τέλος).

Για περισσότερες λεπτομέρειες, ανατρέξτε στις Οδηγίες Χρήσης Magellan.

6.2.4 Εκτέλεση της μεθόδου

Εάν επιλεγεί το Run this method now (Εκτέλεση αυτής της μεθόδου τώρα) στο πλαίσιο διαλόγου Save as (Αποθήκευση ως) στον οδηγό Create/edit a method (Δημιουργία/επεξεργασία μεθόδου), το πλαίσιο διαλόγου Start Measurement Wizard/Start Measurement (Οδηγός έναρξης μέτρησης/Έναρξη μέτρησης) θα εμφανιστεί αφού κάνετε κλικ στο Save (Αποθήκευση). Στο διάλογο Start Measurement (Έναρξη μέτρησης), αντιστοιχίζεται ένα προεπιλεγμένο όνομα χώρου εργασίας, το οποίο ο χρήστης μπορεί να αλλάξει εάν θέλει.

rt Measurement			
Measurement			Measurement parameters
Workspace: Method: Sample ID list: Instrument Use sta Plate in Movements	05112009-001.wsp Method1.mth cker Temp. control Injector control	Arb. cycle kin Modify layout Insert Current: n. def °C Target: n. def °C Optimize Z-position	Plate Plate Description: [GRE96/t] - Greiner 96 Flat Tr. Plate with Cover: No Barcode: No Part of Plate Range: A1:H12 Absorbance Measurement wavelength: 492 nm Reference wavelength: 620 nm Label: Label1
Help Cancel	<<< Back		START
			© 2009 Tecan



Κάντε κλικ στο **Start** (Έναρξη) για να ξεκινήσει η μέτρηση. Θα δημιουργηθεί αυτόματα ένας χώρος εργασίας, ο οποίος περιέχει όλες τις προηγουμένως εισαχθείσες πληροφορίες και ο οποίος θα συλλέξει όλες τις τιμές μέτρησης. Κατά την εκτέλεση της μέτρησης, εμφανίζεται ένα πλαίσιο διαλόγου κατάστασης της μέτρησης που υποδεικνύει την πρόοδο της μέτρησης.

Μετά την ολοκλήρωση της μέτρησης, εμφανίζεται το πλαίσιο διαλόγου **Results** (Αποτελέσματα), στο οποίο προβάλλονται όλα τα αποτελέσματα και οι υπολογισμοί. Υπάρχει η πιθανότητα να παρουσιαστούν σφάλματα κατά την εκτέλεση μέτρησης χωρίς τα αντίστοιχα υγρά (π.χ. πρότυπα).

6.2.5 Αξιολόγηση αποτελέσματος

Επιλέξτε **Evaluate results** (Αξιολόγηση αποτελεσμάτων) για να προβάλετε και αξιολογήσετε τα πρωτογενή δεδομένα. Είναι δυνατή η προβολή των παραμέτρων αξιολόγησης και η επαναξιολόγηση των δεδομένων.

Αυτή η ενότητα σας καθοδηγεί στον οδηγό **Evaluate Results** (Αξιολόγηση αποτελεσμάτων) με τη χρήση ενός αρχείου χώρου εργασίας που εγκαθίσταται αυτόματα με την εγκατάσταση του Magellan.

Στο πλαίσιο διαλόγου *Wizard List (Λίστα οδηγών)*, κάντε κλικ στο **Evaluate** results (Αξιολόγηση αποτελεσμάτων).

Κάντε κλικ στο **Next** (Επόμενο) στη σελίδα **Welcome** (Καλώς ορίσατε) στον οδηγό **Evaluate Results** (Αξιολόγηση αποτελεσμάτων). Εμφανίζεται το πλαίσιο διαλόγου **Select a file** (Επιλογή αρχείου).

Επιλέξτε το χώρο εργασίας Quantitative ELISA

example_Sunrise_InfiniteF50.wsp από τη λίστα και κάντε κλικ στο Make your selection (Επιλογή). Πραγματοποιούνται υπολογισμοί και εμφανίζεται το παρακάτω παράθυρο διάταξης πλάκας:

Evaluate results - Quantitative BLISA example_Surrise_InfiniteF30.wsp - Quantitative BLISA example_Surrise_mth. 🔀															
Fie Ed. Instrument															
	5 1														
10	Ť														
Measurement data 📀 Measurement data Reference data	Evaluate re	0	1	2	3	4	5	6	7	8	9	10	11	12	
Measurement data - Mean	esult		BL1	ST1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37	
v - Measurement data		A	1/1	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Reference data - Mean			0.004	0.207	0.1	0.816	0.174	0.166	0.083	0.085	0.085	80.0	0.162	0.131	
s - Reference data	Edit m	D	NC1	ST1_4	SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
Measurement data as Colors	etho	D	1/2	1/2	0.784	0.212	0.196	0.156	0.106	0.216	0.123	0.103	0.131	0.152	
Reference data as Colors			NC1	ST1 4	SM1 2	SM1 6	SM1 10	SM1_14	SM1 18	SM1 22	SM1 26	SM1 30	SM1 34	SM1 38	
Neduced data 🔗		C	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Difference data			0.068	0.418	0.764	0.205	0.193	0.155	0.104	0.21	0.129	0.112	0.125	0.155	
Difference data - Mean			ST1_1	ST1_5	SM1_3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
v - Difference data			1/2		1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
Difference data as Colors			0.052	0.838	0.64	0.083	0.098	0.108	0.121	0.11	0.162	0.105	0.127	0.093	
Transformed data		e l	ST1_1	ST1_5	SM1_3	SM1_7	SM1_11	SM1_15	SM1_19	SM1_23	SM1_27	SM1_31	SM1_35	SM1_39	
3			2/2	2/2	0.629	2/2	2/2	0.11	0.125	0.115	2/2	0.111	0.120	2/2	
Concentrations S			ST1 2	ST1 6	SM1 4	SM1 8	SM1 12	SM1_16	SM1_20	SM1 24	SM1_28	SM1 32	SM1 36	SM1_40	
Qualitative Results 🛞		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	
🏾 Method layout 🛛 🛞		0	0.103	1.658	0.323	0.104	0.078	0.153	0.143	0.165	0.112	0.094	0.135	0.143	
Of Validation			ST1_2	ST1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	stria
QC Validation criteria		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	Aus
Miscallanaous			0.103	1.655	0.314	0.099	0.079	0.155	0.136	0.164	0.116	0.092	0.124	0.149	CAN
Remarks		P.0	SI1_3	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37		Ë
Error protocol			0 206	0 105	0.845	0 199	0 167	0.085	0.081	0.082	0.081	0 167	0 127		
Audit Trail Signature															
		<u> </u>													
	P	Measurement par	ameters												~
SUMUSE															
		Plate Plate Description:	[SUNRISE_Stan	idard] - Greiner	96 Flat										~
Help	Heb														
Cancel <<< Back														NEXT	
													@ 200	9 Tecan	

Σε κάθε κυψελίδα εμφανίζεται η υπολογισμένη τιμή. Ανάλογα με το επιλεγμένο στοιχείο στη γραμμή ελέγχου, το παράθυρο διάταξης πλάκας αλλάζει αντίστοιχα. Οι παράμετροι και οι ρυθμίσεις μπορούν να αλλαχτούν μέσω των στοιχείων στη γραμμή ελέγχου. Εάν η μέθοδος πρόκειται να τροποποιηθεί, κάντε κλικ στην καρτέλα **Edit method** (Επεξεργασία μεθόδου).



SM1	_10 SM1_14 SM1_18
2	Summary
0.1	Details Edit
SM ²	Edit kinetic settings
1	Copy kinetic settings
0.0	Paste kinetic settings
SM	Graph: Multilabel
2	Graph: Kinetics
0	Graph: Multilabel Kinetics Graph: FLT curves
SM4	Graph: Spectra
	Graph: Dilution series
	Mask/Unmask selection
0.0	Show/Hide layout

Κάντε δεξί κλικ στην κυψελίδα. Εμφανίζεται το σχετικό μενού:

Η επιλογή του στοιχείου **Summary** (Σύνοψη) εμφανίζει το παρακάτω παράθυρο, παρέχοντας αναλυτικές πληροφορίες για τον ορισμό και τις ρυθμίσεις της επιλεγμένης κυψελίδας.

Well: C5	X						
E Method layout							
🗈 📠 Liquid Alias							
🕀 🛲 Replicate							
🕀 📠 Dilution							
🛨 🛲 Multilabel data reduction							
🛨 🛲 Transformation formula(s)							
Kinetic transformation formula(s)							
Concentration transformation formula(s)							
🖭 📠 Strip method names							
🗉 🛄 Instrument data	🗄 🛄 Instrument data						
🗄 💥 Reduced data							
🗄 🔟 Averages							
🕀 🚣 Transformed data							
🗉 🗍 Concentrations							
⊕ 🛕 Statistics							
left Up right							
Expand All	ОК						

Κάντε κλικ στο **Next** (Επόμενο) στο παράθυρο διάταξης πλάκας. Εμφανίζεται το πλαίσιο διαλόγου **Save as** (Αποθήκευση ως), όπου μπορείτε να εισάγετε ένα όνομα αρχείου και σχόλια. Κάντε κλικ στο μικρό κουμπί **Save** (Αποθήκευση) στα αριστερά του παραθύρου ώστε να αποθηκεύσετε το αρχείο. Μπορείτε να συνεχίσετε την εργασία σας στη μέθοδο ή το χώρο εργασίας. Κάντε κλικ στο κουμπί **Finish** (Τέλος) στη δεξιά πλευρά στο κάτω μέρος της οθόνης για να αποθηκεύσετε τον οδηγό. Το πρόγραμμα επιστρέφει στη λίστα οδηγών.



6.2.6 Σύνοψη ορισμού ποσοτικού προσδιορισμού ELISA στο Magellan

1. Αφαίρεση τιμής τυφλών

Ορισμοί στο Magellan

Κάντε κλικ στο Add new transformation (Προσθήκη νέου μετασχηματισμού) στη γραμμή ελέγχου. Εμφανίζεται ένα παράθυρο που σας ρωτά εάν θέλετε να ορίσετε Blank reduction (Αναγωγή απορρόφησης βάσει τυφλών). Κάντε κλικ στο Yes (Ναι) και ο τύπος Blank reduction (Αναγωγή απορρόφησης βάσει τυφλών) αντιστοιχίζεται αυτόματα σε όλες τις κυψελίδες.

2. Ορισμός συγκεντρώσεων

Ορισμοί στο Magellan (Γραμμή ελέγχου – Method layout/ Conc.-, Dil.-, Ref.-values [Διάταξη μεθόδου/τιμές Συγκέντρωσης, Αραίωσης, Αναφοράς])

Επιλεγμένο αναγνωριστικό: ST

Μονάδα: UA/ml

ST1_1	5	(ST1_1Πρότυπο 1, πρώτη πειραματική ομάδα)
ST1_2	10	(ST1_2Πρότυπο 2, πρώτη πειραματική ομάδα)
ST1_3	20	(ST1_3Πρότυπο 3, πρώτη πειραματική ομάδα)
ST1_4	40	(ST1_4Πρότυπο 4, πρώτη πειραματική ομάδα)
ST1_5	80	(ST1_5Πρότυπο 5, πρώτη πειραματική ομάδα)
ST1_6	160	(ST1_6Πρότυπο 6, πρώτη πειραματική ομάδα)

3. Ορισμός πρότυπης καμπύλης

Ορισμοί στο Magellan (Γραμμή ελέγχου – Concentrations/ Standard curve [Συγκεντρώσεις/ Πρότυπη καμπύλη])

Δεδομένα εισόδου αναγωγή απορρόφησης βάσει τυφλών

Τύπος ανάλυσης γραμμική παλινδρόμηση

Άξονας Χ	γραμμικός

Άξονας Υ γραμμικός

4. Ορισμός ορίων

Ορισμοί στο Magellan (Γραμμή ελέγχου – Evaluate data/ Cutoff definition [Αξιολόγηση δεδομένων/ Ορισμός ορίων])

Δεδομένα εισόδου: Μέση συγκέντρωση (UA/ml)

Όρια: 22 18 Θετικό ≥ 22 > ενδιάμεσο ≥ 18 > αρνητικό Μη-ανταγωνιστικός έλεγχος



5. Επικύρωση ελέγχου ποιότητας

Ορισμοί στο Magellan (Γραμμή ελέγχου – Evaluate data/ QC validation [Αξιολόγηση δεδομένων/ Επικύρωση ελέγχου ποιότητας])

Δεδομένα εισόδου: Μία συγκέντρωση (UA/ml)

Συνθήκη επικύρωσης 1 NC1_1<8

Συνθήκη επικύρωσης 2 NC1_2<8

NC1_1.....Αρνητικός μάρτυρας, πρώτο αντίγραφο, πρώτη πειραματική ομάδα

NC1_2.....Αρνητικός μάρτυρας, δεύτερο αντίγραφο, πρώτη πειραματική ομάδα



7. Καθαρισμός, συντήρηση και απόρριψη

7.1 Εισαγωγή

Αυτό το κεφάλαιο περιλαμβάνει τις παρακάτω διαδικασίες:

- Καθαρισμός του οργάνου
- Απολύμανση του οργάνου
- Συντήρηση του οργάνου
- Τοποθέτηση ή αντικατάσταση φίλτρων στον τροχό φίλτρων
- Οδηγίες απόρριψης



ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΠΡΙΝ ΤΟΝ ΚΑΘΑΡΙΣΜΟ Ή ΤΗ ΣΥΝΤΗΡΗΣΗ, ΑΦΑΙΡΕΣΤΕ ΤΗ ΜΙΚΡΟΠΛΑΚΑ.



ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΠΡΙΝ ΤΟΝ ΚΑΘΑΡΙΣΜΟ ΚΑΙ ΤΗΝ ΑΠΟΛΥΜΑΝΣΗ, ΑΠΟΣΥΝΔΕΣΤΕ ΤΟ ΟΡΓΑΝΟ ΑΠΟ ΤΟ ΕΞΩΤΕΡΙΚΟ ΤΡΟΦΟΔΟΤΙΚΟ.



ΠΡΟΣΟΧΗ

ΜΗΝ ΜΕΤΑΚΙΝΕΙΤΕ ΤΟ ΦΟΡΕΑ ΜΕΤΑΦΟΡΑΣ ΤΗΣ ΠΛΑΚΑΣ ΧΕΙΡΟΚΙΝΗΤΑ ΕΚΤΟΣ ΕΑΝ ΤΟ ΟΡΓΑΝΟ ΕΙΝΑΙ ΑΠΕΝΕΡΓΟΠΟΙΗΜΕΝΟ.

7.2 Καθαρισμός του οργάνου



ΠΡΟΕΙΔΟΠΟΙΗΣΗ

Η ΔΙΑΔΙΚΑΣΙΑ ΚΑΘΑΡΙΣΜΟΥ ΠΡΕΠΕΙ ΝΑ ΕΚΤΕΛΕΙΤΑΙ ΣΕ ΑΙΘΟΥΣΑ ΜΕ ΚΑΛΟ ΕΞΑΕΡΙΣΜΟ ΑΠΟ ΕΞΟΥΣΙΟΔΟΤΗΜΕΝΟ ΕΚΠΑΙΔΕΥΜΕΝΟ ΠΡΟΣΩΠΙΚΟ ΠΟΥ ΦΟΡΑΕΙ ΓΑΝΤΙΑ ΜΙΑΣ ΧΡΗΣΗΣ ΚΑΙ ΠΡΟΣΤΑΤΕΥΤΙΚΑ ΓΥΑΛΙΑ ΚΑΙ ΡΟΥΧΙΣΜΟ.

Καθαρίστε το περίβλημα της συσκευής και το φορέα μεταφοράς της πλάκας μόνο με ένα στεγνό ή ελαφρά νοτισμένο ύφασμα. Εάν είναι πολύ βρώμικο, καθαρίστε το με ένα ύφασμα νοτισμένο με έως 70% αιθανόλη ή ήπιο απορρυπαντικό, Microcide SQ ή Decon 90. Στεγνώστε το με χαρτί κουζίνας.

Εάν χυθεί οποιοδήποτε υγρό στο όργανο, θα πρέπει να αφαιρεθεί αμέσως ώστε να αποφευχθεί η διαρροή του υγρού στο σύστημα οπτικών στοιχείων που θα έχει ως αποτέλεσμα απώλεια απόδοσης ή σφάλμα.

7.3 Απολύμανση οργάνου



ΠΡΟΕΙΔΟΠΟΙΗΣΗ

ΕΑΝ ΤΟ ΥΓΡΟ ΠΟΥ ΕΧΕΙ ΧΥΘΕΙ ΣΤΟ ΦΟΡΕΑ ΜΕΤΑΦΟΡΑΣ ΤΗΣ ΠΛΑΚΑΣ ΕΝΔΕΧΕΤΑΙ ΝΑ ΕΙΝΑΙ ΜΟΛΥΣΜΑΤΙΚΟ, ΤΟ ΟΡΓΑΝΟ ΘΑ ΠΡΕΠΕΙ ΝΑ ΑΠΟΛΥΜΑΝΘΕΙ ΣΥΜΦΩΝΑ ΜΕ ΤΟΥΣ ΣΧΕΤΙΚΟΥΣ ΕΘΝΙΚΟΥΣ ΝΟΜΟΥΣ ΚΑΙ ΚΑΝΟΝΙΣΜΟΥΣ.

Όλα τα μέρη του οργάνου που θα έρθουν σε επαφή με βιολογικά δείγματα, δείγματα ασθενών, δείγματα θετικών μαρτύρων ή επικίνδυνο υλικό θα πρέπει να αντιμετωπίζονται ως πιθανόν μολυσματικές περιοχές.



ΠΡΟΕΙΔΟΠΟΙΗΣΗ Η ΔΙΑΔΙΚΑΣΙΑ ΑΠΟΛΥΜΑΝΣΗΣ ΚΑΙ ΤΑ ΑΠΟΛΥΜΑΝΤΙΚΑ ΘΑ ΠΡΕΠΕΙ ΝΑ ΣΥΜΜΟΡΦΩΝΟΝΤΑΙ ΜΕ ΤΟΥΣ ΣΧΕΤΙΚΟΥΣ ΕΘΝΙΚΟΥΣ ΝΟΜΟΥΣ ΚΑΙ ΚΑΝΟΝΙΣΜΟΥΣ.



ΠΡΟΕΙΔΟΠΟΙΗΣΗ ΕΙΝΑΙ ΠΟΛΥ ΣΗΜΑΝΤΙΚΟ ΝΑ ΑΠΟΛΥΜΑΝΘΕΙ ΤΟ ΟΡΓΑΝΟ ΠΛΗΡΩΣ ΠΡΙΝ ΜΕΤΑΦΕΡΘΕΙ ΑΠΟ ΤΟ ΕΡΓΑΣΤΗΡΙΟ Ή ΠΡΙΝ ΠΡΑΓΜΑΤΟΠΟΙΗΘΕΙ ΟΠΟΙΑΔΗΠΟΤΕ ΕΡΓΑΣΙΑ ΕΠΙΣΚΕΥΗΣ/ΣΥΝΤΗΡΗΣΗΣ.

Πριν το όργανο επιστραφεί στον τοπικό αντιπρόσωπο πωλήσεων ή σε κέντρο επισκευής, όλες οι επιφάνειες και ο φορέας μεταφοράς της πλάκας θα πρέπει να απολυμανθούν, ενώ θα πρέπει να συμπληρωθεί και πιστοποιητικό ασφάλειας από το φορέα που το χρησιμοποιεί. Εάν δεν παρασχεθεί πιστοποιητικό ασφάλειας, ο τοπικός αντιπρόσωπος πωλήσεων ή το κέντρο επισκευής ενδέχεται να μην δεχτούν το όργανο ή αυτό να παρακρατηθεί από τις τελωνειακές αρχές.

7.3.1 Διαλύματα απολύμανσης

Οι εξωτερικές επιφάνειες και ο φορέας μεταφοράς της πλάκας θα πρέπει να απολυμανθούν με τη χρήση απολυμαντικών, όπως:

- Microcide SQ
- Decon 90
- Αιθανόλη 70%



ΚΙΝΔΥΝΟΣ ΠΥΡΚΑΓΙΑΣ ΚΑΙ ΕΚΡΗΞΗΣ! ΑΛΚΟΟΛΕΣ ΟΠΩΣ Η ΑΙΘΑΝΟΛΗ ΚΑΙ Η ΙΣΟΠΡΟΠΑΝΟΛΗ ΕΙΝΑΙ ΕΥΦΛΕΚΤΕΣ ΚΑΙ Ο ΛΑΝΘΑΣΜΕΝΟΣ ΧΕΙΡΙΣΜΟΣ ΤΟΥΣ ΜΠΟΡΕΙ ΝΑ ΟΔΗΓΗΣΕΙ ΣΕ ΕΚΡΗΞΕΙΣ Ή/ΚΑΙ ΠΥΡΚΑΓΙΑ. ΠΡΕΠΕΙ ΝΑ ΤΗΡΗΘΟΥΝ ΤΑ ΑΠΑΡΑΙΤΗΤΑ ΜΕΤΡΑ ΑΣΦΑΛΕΙΑΣ ΕΡΓΑΣΤΗΡΙΟΥ.

ΠΡΟΕΙΔΟΠΟΙΗΣΗ



ΠΡΟΣΟΧΗ ΜΗ ΧΡΗΣΙΜΟΠΟΙΕΙΤΕ ΑΚΕΤΟΝΗ ΚΑΘΩΣ ΘΑ ΚΑΤΑΣΤΡΕΨΕΙ ΤΑ ΚΑΛΥΜΜΑΤΑ.



7. Καθαρισμός, συντήρηση και απόρριψη

7.3.2 Διαδικασία απολύμανσης

Εάν το εργαστήριο δεν διαθέτει συγκεκριμένη διαδικασία απολύμανσης, θα πρέπει να χρησιμοποιηθεί η παρακάτω διαδικασία για την απολύμανση των εξωτερικών επιφανειών και του φορέα μεταφοράς της πλάκας του οργάνου.







ΠΡΟΕΙΔΟΠΟΙΗΣΗ Η ΔΙΑΔΙΚΑΣΙΑ ΑΠΟΛΥΜΑΝΣΗΣ ΠΡΕΠΕΙ ΝΑ ΕΚΤΕΛΕΙΤΑΙ ΣΕ ΑΙΘΟΥΣΑ ΜΕ ΚΑΛΟ ΕΞΑΕΡΙΣΜΟ ΑΠΟ ΕΞΟΥΣΙΟΔΟΤΗΜΕΝΟ ΕΚΠΑΙΔΕΥΜΕΝΟ ΠΡΟΣΩΠΙΚΟ ΠΟΥ ΦΟΡΑΕΙ ΓΑΝΤΙΑ ΜΙΑΣ ΧΡΗΣΗΣ ΚΑΙ ΠΡΟΣΤΑΤΕΥΤΙΚΑ ΓΥΑΛΙΑ ΚΑΙ ΡΟΥΧΙΣΜΟ.

ΠΡΟΣΟΧΗ

ΤΟ ΑΠΟΛΥΜΑΝΤΙΚΟ ΓΙΑ ΤΙΣ ΕΠΙΦΑΝΕΙΕΣ ΜΠΟΡΕΙ ΝΑ ΕΠΗΡΕΑΣΕΙ ΑΡΝΗΤΙΚΑ ΤΗΝ ΑΠΟΔΟΣΗ ΤΟΥ ΟΡΓΑΝΟΥ ΕΑΝ ΕΦΑΡΜΟΣΤΕΙ Ή ΕΙΣΕΛΘΕΙ ΚΑΤΑ ΛΑΘΟΣ ΣΤΟ ΟΡΓΑΝΟ.

ΠΡΟΕΙΔΟΠΟΙΗΣΗ

ΠΡΙΝ ΤΗΝ ΑΠΟΛΥΜΑΝΣΗ, ΑΠΟΣΥΝΔΕΣΤΕ ΤΟ ΟΡΓΑΝΟ ΑΠΟ ΤΗΝ ΚΥΡΙΑ ΤΡΟΦΟΔΟΣΙΑ ΩΣΤΕ ΝΑ ΑΠΟΦΥΓΕΤΕ ΤΟΝ ΚΙΝΔΥΝΟ ΠΥΡΚΑΓΙΑΣ Ή ΕΚΡΗΞΗΣ.

- 1. Φοράτε προστατευτικά γάντια, γυαλιά και ρουχισμό.
- Προετοιμάστε ένα κατάλληλο δοχείο για όλα τα υλικά προς απόρριψη που θα χρησιμοποιηθούν κατά τη διαδικασία απολύμανσης.
- 3. Αποσυνδέστε το όργανο από το λογισμικό και απενεργοποιήστε το.
- 4. Αφαιρέστε προσεκτικά το φορέα μεταφοράς της πλάκας.



- 5. Εφαρμόστε προσεκτικά το διάλυμα απολύμανσης στο φορέα μεταφοράς της πλάκας, σύμφωνα με τις Οδηγίες Χρήσης του κατασκευαστή. Μη χρησιμοποιείτε υπερβολική ποσότητα απολυμαντικού ώστε να αποφευχθεί η εισροή του διαλύματος στο όργανο ή το λέρωμα των φακών κατά την τοποθέτηση του φορέα μεταφοράς της πλάκας στη συσκευή.
- 6. Μετά τον απαιτούμενο χρόνο επαφής (σύμφωνα με τις Οδηγίες Χρήσης του κατασκευαστή), σκουπίστε το φορέα μεταφοράς της πλάκας με απαλό χαρτί κουζίνας, νοτισμένο με ένα ήπιο απορρυπαντικό ή αποσταγμένο νερό ώστε να απομακρυνθούν όλα τα ίχνη του απολυμαντικού.



7. Καθαρισμός, συντήρηση και απόρριψη

7. Τοποθετήστε προσεκτικά το φορέα μεταφοράς της πλάκας στο όργανο.



- Εφαρμόστε προσεκτικά το απολυμαντικό στην πλάκα βάσης του φορέα μεταφοράς.
- 9. Μετά τον απαιτούμενο χρόνο επαφής, σκουπίστε την πλάκα βάσης του φορέα μεταφοράς με απαλό χαρτί κουζίνας, νοτισμένο με ένα ήπιο απορρυπαντικό ή αποσταγμένο νερό ώστε να απομακρυνθούν όλα τα ίχνη του απολυμαντικού.
- 10.Εφαρμόστε προσεκτικά το διάλυμα απολυμαντικού σε όλες τις εξωτερικές επιφάνειες του οργάνου.
- 11.Μετά τον απαιτούμενο χρόνο επαφής, σκουπίστε το όργανο με απαλό χαρτί κουζίνας, νοτισμένο με ένα ήπιο απορρυπαντικό ή αποσταγμένο νερό ώστε να απομακρυνθούν όλα τα ίχνη του απολυμαντικού.
- 12. Στεγνώστε την εξωτερική επιφάνεια του οργάνου με απαλό χαρτί κουζίνας.
- 13.Επαναλάβετε τη διαδικασία απολύμανσης σε οποιαδήποτε παρελκόμενα πρόκειται να αφαιρεθούν ή επιστραφούν.
- 14. Απορρίψτε το δοχείο με τα αναλώσιμα σύμφωνα με τους σχετικούς εθνικούς νόμους και κανονισμούς.
- 15.Απολυμάνετε τα χέρια σας και καθαρίστε τα με ένα ήπιο απορρυπαντικό.

Κατά την αποστολή του οργάνου στον τοπικό αντιπρόσωπο πωλήσεων/κέντρο επισκευής, ακολουθήστε τα παρακάτω βήματα:

- 16.Συσκευάστε το όργανο και τα παρελκόμενα.
- 17.Συμπληρώστε το πιστοποιητικό ασφάλειας (δείτε παρακάτω) και προσαρτήστε το πάνω στη συσκευασία έτσι ώστε να είναι εύκολα ορατό.

7.3.3 Πιστοποιητικό ασφάλειας

Για την εξασφάλιση της ασφάλειας και της υγιεινής του προσωπικού, παρακαλούμε τους πελάτες μας να συμπληρώσουν ένα Πιστοποιητικό ασφάλειας (το οποίο παρέχεται με το όργανο) και να τοποθετήσουν ένα αντίγραφο στο πάνω μέρος της συσκευασίας επιστροφής του οργάνου (ορατό στο εξωτερικό μέρος της συσκευασίας!) και ένα δεύτερο αντίγραφο στα έγγραφα αποστολής, πριν την αποστολή στο κέντρο επισκευής για συντήρηση ή επισκευή.

Το όργανο πρέπει να απολυμανθεί στις εγκαταστάσεις του φορέα που το χρησιμοποιεί πριν την αποστολή του (ανατρέξτε στο κεφάλαιο 7.3.2 Ασφάλεια Διαδικασία απολύμανσης).

Η διαδικασία απολύμανσης πρέπει να πραγματοποιηθεί σε μια αίθουσα με καλό εξαερισμό και από εξουσιοδοτημένο και εκπαιδευμένο προσωπικό που θα φοράει γάντια μίας χρήσης χωρίς πούδρα και προστατευτικά γυαλιά και ρουχισμό.

Η διαδικασία απολύμανσης θα πρέπει να πραγματοποιηθεί σύμφωνα με τους εθνικούς, περιφερειακούς και τοπικούς κανονισμούς.



Εάν δεν υποβληθεί πιστοποιητικό ασφάλειας, το κέντρο επισκευής ενδέχεται να μην δεχτεί το όργανο.

Εάν χρειάζεται, το τοπικό τμήμα εξυπηρέτησης πελατών της Tecan μπορεί να σας στείλει ένα νέο αντίγραφο του πιστοποιητικού ασφάλειας.

7.4 Πρόγραμμα προληπτικής συντήρησης για το INFINITE F50 PLUS

Προτείνονται οι παρακάτω διαδικασίες προληπτικής συντήρησης.

7.4.1 Μηνιαία

Καθαρίζετε το περίβλημα και το φορέα μεταφοράς της πλάκας με ένα ήπιο απορρυπαντικό τουλάχιστον μία φορά το μήνα και πιο συχνά εάν κρίνεται απαραίτητο.



ΠΡΟΣΟΧΗ ΜΗ ΧΡΗΣΙΜΟΠΟΙΕΙΤΕ ΑΚΕΤΟΝΗ ΚΑΘΩΣ ΘΑ ΚΑΤΑΣΤΡΕΨΕΙ ΤΑ ΚΑΛΥΜΜΑΤΑ.

7.4.2 Κάθε 4 έτη

Συνιστάται η αντικατάσταση των φίλτρων κάθε 4 έτη.

7.5 Αντικατάσταση και τοποθέτηση φίλτρων

Για να πραγματοποιηθεί τοποθέτηση ή αντικατάσταση φίλτρων με καθοδήγηση από το λογισμικό, το INFINITE F50 PLUS πρέπει να είναι συνδεδεμένο με το Magellan. Εάν κατά τη διαδικασία χαθεί η σύνδεση λόγω ακούσιας διακοπής της σύνδεσης μεταξύ του οργάνου και του υπολογιστή, η λειτουργία του Magellan θα πρέπει να τερματιστεί και το όργανο να απενεργοποιηθεί. Σε αυτήν την περίπτωση, συνεχίστε τη διαδικασία όπως περιγράφεται παρακάτω. Όταν ολοκληρωθεί η διαδικασία πραγματοποιήστε και πάλι τη σύνδεση επανεκκινώντας τη συσκευή και το Magellan και ορίζοντας τα φίλτρα που μόλις τοποθετήθηκαν.



ΠΡΟΣΟΧΗ ΚΑΤΑ ΤΟ ΧΕΙΡΙΣΜΟ ΤΩΝ ΦΙΛΤΡΩΝ, ΠΡΟΣΕΧΕΤΕ ΝΑ ΜΗ ΓΔΑΡΘΟΥΝ Ή ΛΕΡΩΘΟΥΝ ΜΕ ΔΑΧΤΥΛΙΕΣ Ή ΣΚΟΝΗ.



7.5.1 Διαδικασία αλλαγής φίλτρων

Τα φίλτρα του τυπικού τροχού φίλτρων μπορούν να αντικατασταθούν ή να συμπληρωθούν με την ακόλουθη διαδικασία:

- Στο παράθυρο Wizard list (Λίστα οδηγών), κάντε κλικ στο Miscellaneous (Διάφορα).
- 2. Κάντε κλικ στο Instrument control (Έλεγχος οργάνου).
- 3. Κάντε κλικ στο Define filter slides... (Ορισμός slide φίλτρων...).
- Κάντε κλικ στο Filter switching (Αλλαγή φίλτρων) για να ξεκινήσετε τη διαδικασία.
- 5. Αφαιρέστε τη μικροπλάκα από το φορέα μεταφοράς πλάκας!
- Προσεκτικά δώστε κλίση στο όργανο προς τα πίσω έως ότου ακουμπήσει στην πίσω πλευρά με το κάτω μέρος στραμμένο προς το μέρος σας.
- Αφαιρέστε την πλάκα καλύμματος από το κάτω μέρος του οργάνου βγάζοντας τις πορτοκαλί βίδες.



 Αφαιρέστε τον τροχό φίλτρων που είναι στερεωμένος με μαγνήτη, τραβώντας τον προσεκτικά έξω από το όργανο.



9. Τοποθετήστε τον τροχό φίλτρων σε μια καθαρή και επίπεδη επιφάνεια.


7. Καθαρισμός, συντήρηση και απόρριψη

 Κατά την αντικατάσταση ενός φίλτρου, χρησιμοποιήστε το ειδικό εργαλείο για να το αφαιρέσετε από την υποδοχή.

Για το ειδικό εργαλείο και για διαθέσιμα φίλτρα, επικοινωνήστε με τον τοπικό αντιπρόσωπο πωλήσεων της Tecan.



- 11.Ευθυγραμμίστε το ειδικό εργαλείο με την εγκοπή του δακτυλίου ακινητοποίησης. Περιστρέψτε το εργαλείο και αφαιρέστε το δακτύλιο ακινητοποίησης τραβώντας τον έξω από την υποδοχή του φίλτρου.
- 12.Αναποδογυρίστε τον τροχό έτσι ώστε το φίλτρο να γλιστρά έξω από την υποδοχή. Μη χρησιμοποιείτε το ειδικό εργαλείο για να σπρώξετε τα φίλτρα έξω από την υποδοχή, καθώς ενδέχεται να γδαρθούν.
- 13.Το καινούριο φίλτρο πρέπει να τοποθετηθεί στην υποδοχή με το σωστό προσανατολισμό και με προσοχή να μη γδαρθεί ή λερωθεί με δαχτυλιές.





Σημείωση Βεβαιωθείτε ότι το φίλτρο έχει τοποθετηθεί σωστά.



7. Καθαρισμός, συντήρηση και απόρριψη

14.Τοποθετήστε το δακτύλιο ακινητοποίησης στο άκρο του ειδικού εργαλείο και περιστρέψτε τον έτσι ώστε να μην μπορεί να βγει.



- 15.Με τη χρήση του ειδικού εργαλείου, σπρώξτε το δακτύλιο ακινητοποίησης στην υποδοχή του φίλτρου και πιέστε τον σταθερά στη θέση του.
- 16.Περιστρέψτε το εργαλείο έως ότου η εγκοπή στο δακτύλιο ακινητοποίησης ευθυγραμμιστεί με το άκρο του ειδικού εργαλείου. Αφαιρέστε το εργαλείο.



- 17.Επανατοποθετήστε τον τροχό στην υποδοχή του και σπρώξτε τον έως ότου κουμπώσει μαγνητικά.
- 18.Επανατοποθετήστε την πλάκα καλύμματος στο κάτω μέρος του οργάνου με τις τρεις πορτοκαλί βίδες.
- 19. Γυρίστε το όργανο στην όρθια θέση.
- 20.Κάντε κλικ στο **ΟΚ** για να ολοκληρώσετε τη διαδικασία και αρχικοποιήσετε τον τροχό φίλτρων.
- 21.Ορίστε το φίλτρο που μόλις τοποθετήθηκε (για αναλυτική διαδικασία, ανατρέξτε στο επόμενο κεφάλαιο).



7.5.2 Ορισμός φίλτρων

Στο πλαίσιο διαλόγου **Filter Definition** (Ορισμός φίλτρων) αντιστοιχίστε τα κατάλληλα μήκη κύματος στα νέα φίλτρα εισάγοντας τα νέα μήκη κύματος στις αντίστοιχες θέσεις.

Εάν έχει τοποθετηθεί φίλτρο σε νέα θέση, ενεργοποιήστε την κατάλληλη θέση επιλέγοντας το πλαίσιο ελέγχου και εισάγοντας το κατάλληλο μήκος κύματος.

Κάνοντας κλικ στο **Save** (Αποθήκευση), οι ορισμοί των φίλτρων αποθηκεύονται και τα φίλτρα αρχικοποιούνται.

Μετά την αρχικοποίηση των φίλτρων, το όργανο είναι έτοιμο για μετρήσεις.



Σημείωση Προσέξτε να μην μπερδέψετε τις θέσεις των φίλτρων και τα μήκη κύματος καθώς αυτό θα έχει ως αποτέλεσμα λανθασμένα δεδομένα μέτρησης.

7.6 Απόρριψη

7.6.1 Εισαγωγή

Ακολουθήστε τις διαδικασίες εργαστηρίου για την απόρριψη βιολογικά επικίνδυνων υλικών σύμφωνα με εθνικούς και τοπικούς κανονισμούς.

Σε αυτό το κεφάλαιο παρέχονται οδηγίες για τη σύμφωνη με το νόμο απόρριψη αποβλήτων που προκύπτουν από τη χρήση του INFINITE F50 PLUS.



ΠΡΟΣΟΧΗ ΤΗΡΕΙΤΕ ΟΛΟΥΣ ΤΟΥΣ ΚΟΙΝΟΤΙΚΟΥΣ, ΚΡΑΤΙΚΟΥΣ ΚΑΙ ΤΟΠΙΚΟΥΣ ΠΕΡΙΒΑΛΛΟΝΤΙΚΟΥΣ ΚΑΝΟΝΙΣΜΟΥΣ.

7.6.2 Απόρριψη υλικών συσκευασίας

Το υλικό συσκευασίας αποτελείται από ανακυκλώσιμο υλικό. Εάν δεν σκοπεύετε να το κρατήσετε για μελλοντική χρήση, π.χ. για λόγους μεταφοράς και αποθήκευσης, απορρίψτε το υλικό συσκευασίας σύμφωνα με τους τοπικούς κανονισμούς.

7.6.3 Απόρριψη υλικών λειτουργίας

ΠΡΟΕΙΔΟΠΟΙΗΣΗ

ΒΙΟΛΟΓΙΚΟΙ ΚΙΝΔΥΝΟΙ ΕΝΔΕΧΕΤΑΙ ΝΑ ΣΧΕΤΙΖΟΝΤΑΙ ΜΕ ΤΑ ΑΠΟΒΛΗΤΑ (ΜΙΚΡΟΠΛΑΚΑ) ΤΗΣ ΔΙΑΔΙΚΑΣΙΑΣ ΠΟΥ ΕΚΤΕΛΕΙΤΑΙ ΣΤΟ ΜΕΤΡΗΤΗ ΑΠΟΡΡΟΦΗΤΙΚΟΤΗΤΑΣ INFINITE F50 PLUS.



ΧΕΙΡΙΣΤΕΙΤΕ ΤΗ ΧΡΗΣΙΜΟΠΟΙΗΜΕΝΗ ΜΙΚΡΟΠΛΑΚΑ ΚΑΙ ΑΛΛΑ ΑΝΑΛΩΣΙΜΑ, ΚΑΘΩΣ ΚΑΙ ΟΛΕΣ ΤΙΣ ΧΡΗΣΙΜΟΠΟΙΗΜΕΝΕΣ ΟΥΣΙΕΣ ΣΥΜΦΩΝΑ ΜΕ ΤΙΣ ΟΔΗΓΙΕΣ ΓΙΑ ΚΑΛΕΣ ΕΡΓΑΣΤΗΡΙΑΚΕΣ ΠΡΑΚΤΙΚΕΣ.

ΡΩΤΗΣΤΕ ΣΧΕΤΙΚΑ ΜΕ ΤΑ ΚΑΤΑΛΛΗΛΑ ΣΗΜΕΙΑ ΣΥΛΛΟΓΗΣ ΚΑΙ ΤΙΣ ΕΓΚΕΚΡΙΜΕΝΕΣ ΜΕΘΟΔΟΥΣ ΑΠΟΡΡΙΨΗΣ ΣΤΗ ΧΩΡΑ Ή ΣΤΗΝ ΠΕΡΙΟΧΗ ΣΑΣ.



7.6.4 Απόρριψη του οργάνου

Εάν έχετε ερωτήσεις σχετικά με την απόρριψη της συσκευής, επικοινωνήστε με το τοπικό τμήμα εξυπηρέτησης πελατών της Tecan.

Βαθμός ρύπανσης	2 (IEC/EN 61010-1)
Μέθοδος απόρριψης	Μολυσμένα απόβλητα





ΕΞΟΠΛΙΣΜΟΥ (WEEE) ΑΡΝΗΤΙΚΕΣ ΠΕΡΙΒΑΛΛΟΝΤΙΚΕΣ ΕΠΙΠΤΩΣΕΙΣ ΠΟΥ ΣΧΕΤΙΖΟΝΤΑΙ ΜΕ ΤΟ ΧΕΙΡΙΣΜΟ ΑΠΟΒΛΗΤΩΝ.

ΠΡΟΣΟΧΗ ΟΔΗΓΙΑ 2012/19/ΕΕ ΠΕΡΙ ΑΠΟΒΛΗΤΩΝ ΗΛΕΚΤΡΙΚΟΥ ΚΑΙ ΗΛΕΚΤΡΟΝΙΚΟΥ

- ΜΗ ΧΕΙΡΙΖΕΣΤΕ ΗΛΕΚΤΡΙΚΟ ΚΑΙ ΗΛΕΚΤΡΟΝΙΚΟ ΕΞΟΠΛΙΣΜΟ ΩΣ ΟΙΚΙΑΚΑ ΑΠΟΒΛΗΤΑ.
- ΣΥΛΛΕΓΕΤΕ ΞΕΧΩΡΙΣΤΑ ΤΑ ΑΠΟΒΛΗΤΑ ΑΠΟ ΗΛΕΚΤΡΙΚΟ ΚΑΙ ΗΛΕΚΤΡΟΝΙΚΟ ΕΞΟΠΛΙΣΜΟ

ΠΡΟΕΙΔΟΠΟΙΗΣΗ

ΑΝΑΛΟΓΑ ΜΕ ΤΙΣ ΕΦΑΡΜΟΓΕΣ, ΤΜΗΜΑΤΑ ΤΟΥ INFINITE F50 PLUS ΕΝΔΕΧΕΤΑΙ ΝΑ ΕΧΟΥΝ ΕΡΘΕΙ ΣΕ ΕΠΑΦΗ ΜΕ ΒΙΟΛΟΓΙΚΑ ΕΠΙΚΙΝΔΥΝΟ ΥΛΙΚΟ.

• ΧΕΙΡΙΣΤΕΙΤΕ ΑΥΤΟ ΤΟ ΥΛΙΚΟ ΣΥΜΦΩΝΑ ΜΕ ΤΑ ΙΣΧΥΟΝΤΑ ΠΡΟΤΥΠΑ ΚΑΙ ΚΑΝΟΝΙΣΜΟΥΣ ΑΣΦΑΛΕΙΑΣ.

• ΑΠΟΛΥΜΑΝΕΤΕ ΟΛΑ ΤΑ ΤΜΗΜΑΤΑ ΠΡΙΝ ΤΗΝ ΑΠΟΡΡΙΨΗ.

Για περισσότερες πληροφορίες σχετικά με το προϊόν, επικοινωνήστε με:

Tecan Austria GmbH

Untersbergstrasse 1A

A-5082 Grödig/Salzburg

AUSTRIA/EUROPE

T +43 6246 8933 444

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

www.tecan.com



8. Επίλυση προβλημάτων

8.1 Εισαγωγή

Ο εσωτερικός μικροεπεξεργαστής παρακολουθεί και ελέγχει ηλεκτρονικές λειτουργίες, καθώς και μετρήσεις, χειρισμούς και αποτελέσματα. Εάν ο μικροεπεξεργαστής εντοπίσει μια βλάβη ή μια λανθασμένη διαδικασία λειτουργίας, στον υπολογιστή θα εμφανιστεί ένα μήνυμα σφάλματος.

8.1.1 Πίνακας μηνυμάτων σφάλματος και επίλυση προβλημάτων

Ο παρακάτω πίνακας παρέχει μια σύντομη περιγραφή των μηνυμάτων σφάλματος και των ενεργειών επίλυσης προβλημάτων.



Σημείωση Εάν εμφανιστούν μηνύματα σφάλματος που δεν αναφέρονται στον παρακάτω πίνακα, επικοινωνήστε με τον τοπικό αντιπρόσωπο τεχνικής υποστήριξης της Tecan.

Μήνυμα σφάλματος	Περιγραφή	Επίλυση προβλημάτων
Σφάλμα συστήματος		
"Lid Open Error" (Σφάλμα ανοιχτού καλύμματος)	Κάλυμμα ανοιχτό κατά την έναρξη μέτρησης	Κλείστε το κάλυμμα και επανεκκινήστε τη μέτρηση
"MTP Init Error" (Σφάλμα αρχικοποίησης μικροπλάκας))	Δεν ήταν δυνατή η αρχικοποίηση του φορέα μεταφοράς μικροπλάκας	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, σπασμένος ιμάντας ή μηχανικό μπλοκάρισμα φορέα μεταφοράς μικροπλάκας
"MTP lost steps abs(steploss) > max_steploss" (Χαμένα βήματα μικροπλάκας (αρ. βημάτων) > μέγιστος αριθμός)	Χαμένα βήματα μικροπλάκας κατά τη μέτρηση Steploss: αριθμός χαμένων βημάτων max_steploss: αριθμός χαμένων βημάτων που επιτρέπονται	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, δυσκολία μηχανικής κίνησης
"Filter lost steps abs(steploss) > max_steploss" (Χαμένα βήματα φίλτρου (αρ. βημάτων) > μέγιστος αριθμός)	Ο τροχός φίλτρων έχασε βήματα κατά τη μέτρηση. Steploss: αριθμός χαμένων βημάτων max_steploss: αριθμός χαμένων βημάτων που επιτρέπονται	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, δυσκολία μηχανικής κίνησης
"USB timeout" (Λήξη χρόνου αναμονής USB)	Λήξη χρόνου αναμονής στην επικοινωνία με USB	Σφάλμα συστήματος – αναφέρετε στο τμήμα εξυπηρέτησης πελατών
"Lamp Low! Minimum: <i>minimum</i> , Maximum: <i>maximum</i> " (Χαμηλή απόδοση λυχνίας! Ελάχιστο: ελάχιστο, Μέγιστο: μέγιστο) Αυτό το μήνυμα εμφανίζεται με εκδόσεις υλικολογισμικού έως την V1.11.	Η ένταση φωτός που μετρήθηκε δεν έφτασε την αναμενόμενη τιμή μεταξύ ελάχιστου και μέγιστου	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, σπασμένη ίνα



8. Επίλυση προβλημάτων

Μήνυμα σφάλματος	Περιγραφή	Επίλυση προβλημάτων
Σφάλμα συστήματος		
Από την έκδοση υλικολογισμικού V1.12 και ύστερα, το μήνυμα "Lamp Low" (Χαμηλή απόδοση λυχνίας) αντικαθίσταται από το: "Prepare REF check" (Προετοιμασία ελέγχου τιμών αναφοράς)	Η ένταση φωτός που μετρήθηκε δεν έφτασε την αναμενόμενη τιμή μεταξύ <i>minimum</i> (ελάχιστου) και <i>maximum</i> (μέγιστου)	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, σπασμένη ίνα
"Wavelength Not Available ! Wavelength: <i>wavelength</i> nm" (Μήκος κύματος μη διαθέσιμο! Μήκος κύματος: μήκος κύματος nm)	Το φίλτρο με το αναφερόμενο μήκος κύματος δεν βρέθηκε στον τροχό φίλτρων	Σφάλμα συστήματος – αναφέρετε στο τμήμα εξυπηρέτησης πελατών
"Channel Low! Channel: channel_nr, Minimum: minimum, Maximum maximum (Σήμα καναλιού χαμηλό! Κανάλι: αρ. καναλιού, Ελάχιστο: ελάχιστο, Μέγιστο: μέγιστο)	Το σήμα στο κανάλι που αναφέρεται δεν έλαβε την αναμενόμενη τιμή μεταξύ ελάχιστου και μέγιστου	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, σπασμένη ίνα
"Invalid Wavelength! Wavelength: <i>wavelength</i> nm" (Μη έγκυρο μήκος κύματος! Μήκος κύματος: μήκος κύματος nm)	Το μήκος κύματος φίλτρου είναι εκτός του εύρους μηκών κύματος του λευκού και του μπλε LED	Σφάλμα συστήματος – αναφέρετε στο τμήμα εξυπηρέτησης πελατών
"Lamp Overflow!" Minimum: minimum, Maximum: maximum" (Υπερχείλιση λυχνίας! Ελάχιστο: ελάχιστο, Μέγιστο: μέγιστο)	Το σήμα στο μετατροπέα αναλογικού/ψηφιακού (ADC) υπερβαίνει την αναμενόμενη περιοχή μεταξύ ελάχιστου και μέγιστου	Πρόβλημα υλικού: Ηλεκτρονική βλάβη
"Value Not Set: <i>value -1</i> " (Δεν έχει οριστεί τιμή: τιμή -1)	Η αναφερόμενη τιμή δεν έχει οριστεί	Σφάλμα συστήματος – αναφέρετε στο τμήμα εξυπηρέτησης πελατών
"Filter Init Error" (Σφάλμα αρχικοποίησης φίλτρου)	Δεν ήταν δυνατή η αρχικοποίηση του φορέα μεταφοράς φίλτρου	Πρόβλημα υλικού: Ηλεκτρονική βλάβη, μηχανικό μπλοκάρισμα φορέα μεταφοράς τροχού φίλτρων

8.1.2 Ορισμός «υπερχείλισης»

Εάν το αποτέλεσμα μιας μέτρησης απορρόφησης δεν είναι εντός των προδιαγραφών του οργάνου (> 4,0 OD), θα υπάρξει υπερχείλιση και η μετρημένη τιμή OD της αντίστοιχης κυψελίδας θα αντικατασταθεί με το 'Overflow' (Υπερχείλιση). Αυτό γίνεται από το λογισμικό και όχι από το όργανο.

8.1.3 Διακοπή τροφοδοσίας

Σε περίπτωση διακοπής τροφοδοσίας συμβαίνουν τα παρακάτω:

Διακοπή τροφοδοσίας στο όργανο και όχι στον υπολογιστή (π.χ. υπολογιστής συνδεδεμένος σε σύστημα αδιάλειπτης παροχής τροφοδοσίας): η σύνδεση USB μεταξύ του οργάνου και υπολογιστή χάνεται. Δημιουργείται μήνυμα σφάλματος από το λογισμικό.

Διακοπή τροφοδοσίας τόσο στο όργανο όσο και στον υπολογιστή: πρέπει να γίνει επανεκκίνηση του υπολογιστή. Δεδομένα μέτρησης δεν θα είναι διαθέσιμα.



Συντομογραφίες, εμπορικά σήματα και σύμβολα

9.1 Συντομογραφίες

Οι παρακάτω συντομογραφίες παρέχονται ως αναφορά και ενδέχεται να εμφανίζονται στις Οδηγίες Χρήσης.

А	Ampere (Αμπέρ)
AC	Alternating Current (Εναλλασσόμενο ρεύμα)
ADC	Analog Digital Converter (Μετατροπέας αναλογικού/ψηφιακού)
ANSI/SBS	American National Standards Institute/Society for Biomolecular Screening (Αμερικανικό Εθνικό Ινστιτούτο Προτύπων/Εταιρεία Βιομοριακού Ελέγχου)
ASCII	American Standard Code for Information Interchange (Αμερικανικός Πρότυπος Κώδικας για Ανταλλαγή Πληροφοριών)
ASTM	American Society for Testing and Material (Αμερικανικός Οργανισμός Δοκιμών και Υλικών)
°C	Βαθμοί Κελσίου
CE	CE conformity marking – Σήμανση συμμόρφωσης CE
CFR	Code of Federal Regulations (Κώδικας Ομοσπονδιακών Κανονισμών)
cm	Εκατοστά
DC	Direct Current (Συνεχές ρεύμα)
EC	European Community (Ευρωπαϊκή Ένωση)
ELISA	Enzyme-linked Immunosorbent Assay (Ενζυμική ανοσοπροσροφητική μέθοδος προσδιορισμού)
EN	European Norm (Ευρωπαϊκή Νόρμα)
°F	Βαθμοί Φαρενάιτ
FDA	Food and Drug Administration (Οργανισμός Τροφίμων και Φαρμάκων)
Hz	Hertz
IEC	International Electrotechnical Commission (Διεθνής Ηλεκτροτεχνική Επιτροπή)
ID	Identification (Αναγνωριστικό)
IFU	Instructions for Use (Οδηγίες Χρήσης)
IQ	Installation Qualification (Επαλήθευση εγκατάστασης)
IVD	In vitro diagnostics (Διαγνωστικός έλεγχος in vitro)
IVDR	In Vitro Diagnostic Regulation (IVDR) (EU) 2017/746 Κανονισμός In Vitro Διάγνωσης (IVDR) (EE) 2017/746
kg	Κιλά
1	Λίτρα
LED	Light Emitting Diode (Δίοδος εκπομπής φωτός)
LIS	Laboratory Information System (Πληροφοριακό Σύστημα Εργαστηρίου)
mg	Χιλιοστογραμμάρια
ml	Χιλιοστόλιτρα



9. Συντομογραφίες, εμπορικά σήματα και σύμβολα

mm	Χιλιοστά
MTP	Microplate (Μικροπλάκα)
μΙ	Μικρολίτρα
NFM	Nonferrous Metal (Μη σιδηρούχα μέταλλα)
NIST	National Institute of Standards and Technology (Εθνικό Ινστιτούτο Προτύπων και Τεχνολογίας)
nm	Νανόμετρα
NRTL	Nationally Recognized Testing Laboratory (Εθνικά Αναγνωρισμένο Εργαστήριο Δοκιμών)
OD	Optical Density (Οπτική πυκνότητα)
OQ	Operational Qualification (Επαλήθευση λειτουργικότητας)
PCB	Printed Circuit Board (Διοικητικό Συμβούλιο τυπωμένων κυκλωμάτων)
RF	Radio Frequency (Ραδιοσυχνότητα)
RoHS	Restriction of the Use of Certain Hazardous Substances (Περιορισμός χρήσης ορισμένων επικίνδυνων ουσιών)
SOP	Standard Operating Procedure (Πρότυπη διαδικασία λειτουργίας)
USB	Universal Serial Bus (Ενιαίος Σειριακός Δίαυλος)
UA	Arbitrary Units (Σχετικές μονάδες)
TÜV	Technischer άberwachungsverein (Οργανισμός τεχνικών επιθεωρήσεων)
V	Volt (Βολτ)
VA	Volt Ampere (Βολταμπέρ)
WEEE	Waste electrical and electronic equipment (Απόβλητα ηλεκτρικού και ηλεκτρονικού εξοπλισμού)

9.2 Εμπορικά σήματα

Τα παρακάτω ονόματα προϊόντων και οποιαδήποτε κατατεθέντα ή μη εμπορικά σήματα που αναφέρονται στο παρόν έγγραφο χρησιμοποιούνται μόνο για σκοπούς ταυτοποίησης και αποτελούν την αποκλειστική ιδιοκτησία των αντίστοιχων κατόχων τους.

- Οι επωνυμίες Magellan[™], Infinite[®], MultiCheck[™], Tecan[®] και το λογότυπο Tecan είναι εμπορικά σήματα της Tecan Group Ltd., Männedorf, Switzerland
- Οι επωνυμίες Windows® και Excel® είναι σήματα κατατεθέντα της Microsoft Corporation, Redmond, WA, USA
- Οι επωνυμίες Pentium[®] και Atom[™] είναι εμπορικά σήματα της Intel Corporation, Santa Clara, CA, USA
- Η επωνυμία Adobe® Reader® είναι σήμα κατατεθέν της Adobe Systems Incorporated, Seattle, WA, USA
- Η επωνυμία Microcide SQ[™] είναι εμπορικό σήμα της Global Biotechnologies Inc., Portland, ME, USA
- Η επωνυμία Decon 90[™] είναι εμπορικό σήμα της Decon Laboratories Ltd., Hove, East Sussex, UK



9.3 Σύμβολα

Τα παρακάτω σύμβολα εμφανίζονται στο όργανο.

	Κατασκευαστής
[]	Ημερομηνία κατασκευής
(€	Σήμανση CE
UK CA	Αξιολόγηση της συμμόρφωσης Ηνωμένου Βασιλείου Η παρούσα σήμανση δείχνει ότι το προϊόν με αυτή την ετικέτα συμμορφώνεται με τον ισχύοντα κανονισμό στη Μεγάλη Βρετανία.
ĺ	Διαβάστε τις Οδηγίες Χρήσης πριν χρησιμοποιήσετε το όργανο
IVD	Ιατροτεχνολογικά βοηθήματα που χρησιμοποιούνται στη διάγνωση in vitro
UDI	Αποκλειστικό αναγνωριστικό τεχνολογικού προϊόντος Το σύμβολο UDI προσδιορίζει τον φορέα δεδομένων στην ετικέτα.
REF	Αριθμός παραγγελίας
SN	Αυξάνων αριθμός
•	Ετικέτα USB
	Σύμβολο WEEE
(50)	Σύμβολο RoHS, Κίνα
	NRTL TÜV SÜD MARK
	Βιολογικά επικίνδυνο



Ευρετήριο

Α

Absorbance (Απορρόφηση)
С
Comment (Σχόλιο)25
I
Incubation (Επώαση)26
К
Kinetic (Ανάλυση κινητικής).
Γραμμή ελέγχου24
Kinetic Condition
(Συνθήκη ανάλυσης κινητικής)
L
Lab Ware (Εργαστήριο), Γραμμή ελέγχου 22
М
Magellan Tracker 17
Measurements (Μετρήσεις),
Γραμμή ελέγχου
Miscellaneous (Διαφορα), Ι ραμμη ελεγχου 25
Ρ
Part of Plate (Τμήμα πλάκας)22 Plate (Πλάκα)22
S
Shaking (Ανακίνηση)
U
User Request (Aímua vonom) 25
W
vvait (Timer) (Αναμονή [Χρονοοιακοπτής]) 25
A
Αλλαγή φίλτρων72
Ανακίνηση35
Άνοιγμα συσκευασίας
Άνοιγμα συσκευασίας και επιθεώρηση13
Διαδικασία ανοίγματος συσκευασίας
Αντικατάσταση φίλτρων
Απαιτήσεις συστήματος15
Απαιτησεις τροφοοοσιας
Αποοοση, βελτιστοποιηση
Απολυμανση
Διαοικασια
Διαλυματα
Γιιστοποιητικό ασφαλειας
- ΤΤΟΡΡΙΨΙ]
Ομγανο

Υλικό συσκευασίας	75
Ασφάλεια	7

В

Βελτιστοποίηση απόδο	σης 33
----------------------	--------

Г

Γραμμή ελέγχου	22
Actions (Ενέργειες)	23
Kinetic (Ανάλυση κινητικής)	24
Lab Ware (Εργαστήριο)	22
Measurements (Μετρήσεις)	23
Miscellaneous (Διάφορα)	25

Δ

Διαδικασία αυτόματου ελέγχου	41
Διακοπή τροφοδοσίας	78

Ε

Έλεγχος ποιότητας	41
Ενεργοποίηση του οργάνου	16
Επαλήθευση λειτουργικότητας	41
Έλεγχος MultiCheck	41
Έλεγχος ακρίβειας	42
Έλεγχος γραμμικότητας	43
Έλεγχος μικροπλάκας	41
Επίλυση προβλημάτων	77

Κ

7
1

Λ

Λίστα οδηγών	19
Λογισμικό	17
Εγκατάσταση	17
Επαλήθευση εγκατάστασης	18
Λίστα οδηγών	19
Περιβάλλον χρήσης	19

М

Μετρήσεις ανάλυσης κινητικής	30
Μετρήσεις πολλαπλής σήμανσης	29
Μετρήσεις τελικού σημείου	28
Μηνύματα σφαλμάτων	77
Μικροπλάκες	39
Βελτιστοποίηση απόδοσης	. 34

0

Οδηγός	
Αξιολόγηση αποτελεσμάτων	
Δημιουργία/επεξεργασία λίστας	
αναγνωριστικών δειγμάτων	
Δημιουργία/επεξεργασία μεθόδου	
Έναρξη μέτρησης	19
Επισύναψη υπογραφής	
Λίστα οδηγών	19



Ευρετήριο

Όργανο	
Θέση	33
Λειτουργίες	35
Παρελκόμενα	40
Περιγραφή	35
Προδιαγραφές	
Ορισμός μετρήσεων	
Ορισμός φίλτρων	75

Π

Παράδειγμα εφαρμογής	45
Ποσοτικός προσδιορισμός ELISA	45
Παράθυρο πληροφοριών	27
Παράθυρο ροής εργασιών	26
Περιβαλλοντικές απαιτήσεις	14
Πιστοποιητικό ασφάλειας	70
Πρόγραμμα επεξεργασίας παραμέτρων	
μέτρησης	21
Προδιαγραφές	38
Προφίλ χρηστών	11

Σ

Στοιχεία προγράμματος	
Absorbance (Απορρόφηση)	23

Comment (Σχόλιο)	25
Incubation (Επώαση)	26
Kinetic Condition (Συνθήκη ανάλυσης κινητικής).	24
Kinetic Cycle (Κύκλος ανάλυσης κινητικής)	24
Part of Plate (Τμήμα πλάκας)	22
Plate (Πλάκα)	22
Shaking (Ανακίνηση)	23
User Request (Αίτημα χρήστη)	25
Wait (Timer) (Αναμονή [Χρονοδιακόπτης])	25
Δέσμευση και αποδέσμευση	31
Σύμβολα	81
Συντήρηση67,	71
Πρόγραμμα συντήρησης	71
Συντομογραφίες	79
Τ	
Τροχός φίλτρων	37
Y	
Υπερχείλιση	78



Εξυπηρέτηση πελατών της Tecan

Εάν έχετε ερωτήσεις ή χρειάζεστε τεχνική υποστήριξη για το προϊόν Tecan που διαθέτετε, επικοινωνήστε με τον τοπικό εκπρόσωπο Εξυπηρέτησης Πελατών της Tecan. Για στοιχεία επικοινωνίας, μεταβείτε στον ιστότοπο <u>http://www.tecan.com/</u>.

Πριν επικοινωνήσετε με την Tecan για την παροχή υποστήριξης για το προϊόν, φροντίστε να έχετε διαθέσιμες τις ακόλουθες πληροφορίες, ώστε να λάβετε την καλύτερη δυνατή τεχνική υποστήριξη (βλ. πλακέτα στοιχείων προϊόντος):

- Όνομα μοντέλου του προϊόντος σας
- Αριθμός σειράς (SN) του προϊόντος σας
- Λογισμικό και έκδοση λογισμικού (αν υπάρχουν)
- Περιγραφή του προβλήματος και στοιχεία υπεύθυνου επικοινωνίας
- Ημερομηνία και ώρα όπου παρουσιάστηκε το πρόβλημα
- Ενέργειες στις οποίες έχετε ήδη προβεί για να διορθώσετε το πρόβλημα
- Τα στοιχεία επικοινωνίας σας (αριθμός τηλεφώνου, αριθμός φαξ, διεύθυνση email, κ.λπ.)

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.