# **•**TECAN•

Руководство по эксплуатации

фотометра

# **INFINITE F50 PLUS**

Программное обеспечение

# MAGELLAN



№ документа:	30186919
	2023-06
Редакция документа:	1.2
Версия микропрограмма:	3.33
Версия ПО <b>Magellan</b> :	7.5







#### ПРЕДУПРЕЖДЕНИЕ

#### ПЕРЕД НАЧАЛОМ ЭКСПЛУАТАЦИИ ОБОРУДОВАНИЯ ВНИМАТЕЛЬНО ПРОЧИТАЙТЕ И ВЫПОЛНИТЕ УКАЗАНИЯ, ПРИВЕДЕННЫЕ В НАСТОЯЩЕМ РУКОВОДСТВЕ.

#### Примечание

При составлении настоящего руководства были приложены все усилия, чтобы избежать ошибок в тексте и на диаграммах, однако компания Tecan Austria GmbH не несет ответственности за какие-либо ошибки, которые могут содержаться в настоящем руководстве.

Стратегия Tecan Austria GmbH предусматривает совершенствование изделий по мере появления новых методик и компонентов. Поэтому Tecan Austria GmbH оставляет за собой право изменения технических характеристик в любое время *при условии выполнения надлежащих требований контроля, проверки и аттестации.* 

Мы будем рады любым комментариям в отношении настоящего документа.



#### Производитель

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

#### Информация об авторском праве

Информация, содержащаяся в настоящем документе, является собственностью компании Tecan Austria GmbH, и ее копирование, тиражирование или передача другому лицу или лицам не допускаются без предварительного письменного разрешения.

Авторское право © Tecan Austria GmbH Все права защищены. Напечатано в Австрии.

#### Заявление о соответствии стандартам ЕС

См. последнюю страницу этого руководства по эксплуатации.

#### О руководстве по эксплуатации

Оригинальное руководство по эксплуатации. Настоящий документ является **Руководством по эксплуатации** (далее — «Руководство») фотометра INFINITE F50 PLUS, предназначенного для измерения поглощения (оптической плотности) образцов на 96-луночных микропланшетах. Руководство содержит справочную информацию и инструкции для пользователя.

В Руководстве содержатся следующие инструкции:

- установка прибора;
- работа с прибором;
- чистка и техническое обслуживание прибора.



#### Замечание к снимкам экранов

Номер версии, показанный на снимках экранов, может не совпадать с номером текущей версии программного обеспечения. Замена снимков экрана производится только при изменении содержания, относящегося к применению.

#### История изменений Руководства по эксплуатации

Версия	Изменения	
V 1.2	Новая DoC (Декларация о соответствии)	

## Предупреждения, предостережения и примечания

В настоящем руководстве используются следующие типы примечаний, подчеркивающих важную информацию или предупреждающих пользователя о потенциально опасной ситуации:

> Примечание Содержит полезные сведения.

ПРЕДОСТЕРЕЖЕНИЕ УКАЗЫВАЕТ НА ВОЗМОЖНОСТЬ ПОВРЕЖДЕНИЯ ПРИБОРА ИЛИ ПОТЕРИ ДАННЫХ В СЛУЧАЕ НЕСОБЛЮДЕНИЯ ИНСТРУКЦИЙ.



**STOP** 





# ОБОРУДОВАНИЯ В СЛУЧАЕ НЕСОБЛЮДЕНИЯ ИНСТРУКЦИЙ.

ПРЕДУПРЕЖДЕНИЕ УКАЗЫВАЕТ НА ВОЗМОЖНОСТЬ ПОЛУЧЕНИЯ ТЯЖЕЛЫХ ТРАВМ, НАСТУПЛЕНИЯ СМЕРТИ ИЛИ ПОВРЕЖДЕНИЯ

ПРЕДУПРЕЖДЕНИЕ

УКАЗЫВАЕТ НА ВОЗМОЖНОЕ НАЛИЧИЕ БИОЛОГИЧЕСКИ ОПАСНОГО МАТЕРИАЛА. В ЛАБОРАТОРИИ НЕОБХОДИМО СОБЛЮДАТЬ СООТВЕТСТВУЮЩИЕ МЕРЫ ТЕХНИКИ БЕЗОПАСНОСТИ.

#### ПРЕДУПРЕЖДЕНИЕ

УКАЗЫВАЕТ НА ВОЗМОЖНОЕ НАЛИЧИЕ ЛЕГКОВОСПЛАМЕНИМЫХ МАТЕРИАЛОВ И ОПАСНОСТЬ ПОЖАРА. В ЛАБОРАТОРИИ НЕОБХОДИМО СОБЛЮДАТЬ СООТВЕТСТВУЮЩИЕ МЕРЫ ТЕХНИКИ БЕЗОПАСНОСТИ.





#### ВНИМАНИЕ

ОТРИЦАТЕЛЬНОЕ ВЛИЯНИЕ НА ОКРУЖАЮЩУЮ СРЕДУ, СВЯЗАННОЕ С ДИРЕКТИВОЙ 2012/19/EU ПО ОБРАБОТКЕ ОТХОДОВ ЭЛЕКТРИЧЕСКОГО И ЭЛЕКТРОННОГО ОБОРУДОВАНИЯ



- (DIRECTIVE 2012/19/EU ON WASTE ELECTRICAL AND ELECTRONIC EQUIPMENT (WEEE)).
- НЕ ВЫБРАСЫВАЙТЕ ЭЛЕКТРИЧЕСКОЕ И ЭЛЕКТРОННОЕ ОБОРУДОВАНИЕ ВМЕСТЕ С НЕСОРТИРОВАННЫМ БЫТОВЫМ МУСОРОМ.
- СОБИРАЙТЕ ОТХОДЫ ЭЛЕКТРИЧЕСКОГО И ЭЛЕКТРОННОГО ОБОРУДОВАНИЯ ОТДЕЛЬНО.



## Содержание

1.	Техн	ника без	зопасности	7
	1.1	Техник	ка безопасности при работе с прибором	7
2.	Обш	цие свед	дения	9
	2.1	Назнач	Чение	
	2.2	Принц	ИП ДЕЙСТВИЯ	
	2.3	Профи	ль пользователя	
		2.3.1	Профессиональный пользователь - уровень администратора	
		2.3.2	Конечный или рядовой пользователь	
		2.3.3	Специалисты по ремонту	
3	При		клаботе	13
<b>J</b> .	2.1	Васпан		
	3.1			
		3.1.1	Осмотр содержимого поставки	دا 12
	2 2	J.I.Z	Порядок распаковки	13 14
	3.Z	Требон	вания к питанию	
	3.3	Требон	вания к окружающей среде Вания к окстомо	
	3.4 2 E	Преоог	вания к системе	
	3.5	Програ	ение приоора	
	3.0	261	Врадониа / Назначание	
		3.0.1	Введение / Пазначение	17
		3.0.2	Порядок установки	
		3.0.3	Проверка установки	
	27	Magall	Пачало расоты с программным осеспечением маденал	
	3.7	371	ан – гедактор параметров измерения	ຊາ ວາ
		372	Панель управления	
		373	Информационная пачель	
	3.8	0.7.0	ипформационная папель	27 27
	5.0	3 8 1	Определение измерении Определение измерений конеценти тоцки	
		382	Определение измерений с несколькими метками	27 28
		383	Определение измерении с несколькими метками	20 20
		384	Создание отступов и разблокирование программных элементов	
	39	0.0.4 Оптим		
	0.0	391	Расположение прибора	
		392	Процедура эксплуатации	
	0	0.0.2	прододура околучации	ວາ
4.	Опи	сание п	риоора	
	4.1	Описа	ние прибора	
		4.1.1	Встряхивание микропланшета	
	4.2	Описа	ние приоора	
	4.3	Описа	ние диска фильтров	
	4.4	Технич	ческие характеристики приоора	
		4.4.1	Оощие характеристики	
		4.4.Z	Характеристики измерения	
	4 E	4.4.3		
-	4.5	прина	длежности приоора	
5.	Кон	гроль ка	ачества	39
	5.1	Введе	ние	39
	5.2	Проце,	дура самопроверки	39
	5.3	Эксплу	уатационная квалификация (ЭК)	39
		5.3.1	Тест MultiCheck	39
		5.3.2	Тестирование микропланшета	39
		5.3.3	Тест сходимости прибора на образцах жидкости	40
		5.3.4	Iест линейности прибора на образцах жидкости	41
6.	При	мер при	именения	43
	6.1	Введе	ние	43
	62	Приме	р с пошаговыми инструкциями: количественный анализ ELISA	43



	6.2.1	Описание диагностического комплекта	43
	6.2.2	Создание метода	45
	6.2.3	Настройка печатного отчета	56
	6.2.4	Запуск метода	60
	6.2.5	Оценка результатов	61
	6.2.6	Определение количественного анализа ELISA	63
7. Чи	стка, техн	ическое обслуживание и утилизация	65
7.1	Введен	ие	65
7.2	Чистка	прибора	65
7.3	Дезинф	рекция прибора	66
	7.3.1	Растворы для дезинфекции	66
	7.3.2	Процедура дезинфекции	67
	7.3.3	Сертификат безопасности	69
7.4	План п	рофилактического технического обслуживания фотометра INFINITE F50 PLU	JS70
	7.4.1	Ежемесячно	70
	7.4.2	Каждые четыре года	70
7.5	Замена	а и установка фильтров	70
	7.5.1	Замена фильтров	71
	7.5.2	Определение фильтров	74
7.6	Утилиз	ация	74
	7.6.1	Введение	74
	7.6.2	Утилизация упаковочных материалов	74
	7.6.3	Утилизация рабочих материалов	74
	7.6.4	Утилизация прибора	75
8. По	иск неисг	<b>травностей</b>	77
8.1	Введен	ие	77
	8.1.1	Таблица сообщений об ошибках и поиск неисправностей	77
	8.1.2	Переполнение	78
	8.1.3	Сбой электропитания	79
9. Co	кращения	а, товарные знаки и символы	81
9.1	Сокрац	цения	81
9.2	Торгов	ые знаки	82
9.3	Симво	ЛЫ	83
Указател	1Ь		85
Сервисн	ые центр	ы Тесап	87



## 1. Техника безопасности

## 1.1 Техника безопасности при работе с прибором

- 1. Для снижения риска травмы, пожара или поражения электрическим током при эксплуатации данного прибора всегда соблюдайте общие правила техники безопасности.
- 2. Внимательно прочтите всю информацию, приведенную в Руководстве. Пренебрежение инструкциями, а также их непонимание или несоблюдение может привести к повреждению прибора, травмированию обслуживающего персонала или неправильной работе прибора. Компания Тесап не несет ответственности за повреждения или травмы, возникшие в результате неправильного обращения с прибором.
- 3. Соблюдайте все инструкции, обозначенные в Руководстве пиктограммами «ПРЕДУПРЕЖДЕНИЕ» и «ПРЕДОСТЕРЕЖЕНИЕ».
- 4. Всегда отключайте прибор от электросети перед техническим обслуживанием, чисткой и дезинфекцией.
- 5. Никогда не вскрывайте корпус прибора.
- Соблюдайте надлежащие правила техники безопасности для лабораторий, такие как ношение защитной спецодежды (перчатки, лабораторный халат, защитные очки) и применение надлежащих одобренных процедур техники безопасности в лабораториях.



#### ПРЕДОСТЕРЕЖЕНИЕ

#### НЕСОБЛЮДЕНИЕ ПРИВЕДЕННЫХ В РУКОВОДСТВЕ ИНСТРУКЦИЙ МОЖЕТ ПРИВЕСТИ К ПОВРЕЖДЕНИЮ ПРИБОРА, НЕШТАТНОМУ РЕЖИМУ ВЫПОЛНЕНИЯ ОПЕРАЦИЙ И НАРУШЕНИЮ БЕЗОПАСНОСТИ ПРИБОРА.

Предполагается, что операторы прибора обладают необходимым профессиональным опытом и хорошо ознакомлены с правилами техники безопасности при обращении с химикатами и веществами, представляющими биологическую опасность.

Соблюдайте следующие законы и рекомендации:

- государственные законы о безопасности в промышленности;
- нормативные правила по предупреждению производственного травматизма;
- инструкции производителей реагентов, содержащиеся в паспортах безопасности этих веществ.







В ЗАВИСИМОСТИ ОТ ПРИМЕНЕНИЯ ПРИБОРА INFINITE F50 PLUS, НА НЕГО МОЖЕТ ПОПАСТЬ ВЕЩЕСТВО, ПРЕДСТАВЛЯЮЩЕЕ ИНФЕКЦИОННУЮ ИЛИ ИНУЮ БИОЛОГИЧЕСКУЮ ОПАСНОСТЬ.

К РАБОТЕ С ПРИБОРОМ ДОЛЖНЫ ДОПУСКАТЬСЯ ТОЛЬКО КВАЛИФИЦИРОВАННЫЕ СПЕЦИАЛИСТЫ. ПЕРЕД ВЫПОЛНЕНИЕМ ЕГО ОБСЛУЖИВАНИЯ, ПЕРЕМЕЩЕНИЯ ИЛИ УТИЛИЗАЦИИ, ВСЕГДА ДЕЗИНФИЦИРУЙТЕ ПРИБОР В СООТВЕТСТВИИ С ИНСТРУКЦИЯМИ, ПРИВЕДЕННЫМИ В НАСТОЯЩЕМ РУКОВОДСТВЕ.

ПРИ ВЫПОЛНЕНИИ РАБОТ С ВЕЩЕСТВАМИ, ПРЕДСТАВЛЯЮЩИМИ ИНФЕКЦИОННУЮ ОПАСНОСТЬ, СОБЛЮДАЙТЕ НАДЛЕЖАЩИЕ ПРАВИЛА ТЕХНИКИ БЕЗОПАСНОСТИ ДЛЯ ЛАБОРАТОРИЙ, ТАКИЕ КАК НОШЕНИЕ ЗАЩИТНОЙ СПЕЦОДЕЖДЫ.

#### ПРЕДУПРЕЖДЕНИЕ

ДАННЫЙ ПРИБОР ОТВЕЧАЕТ ТРЕБОВАНИЯМ ПО ЭМИССИИ И ПОМЕХОУСТОЙЧИВОСТИ, СОДЕРЖАЩИМСЯ В СТАНДАРТЕ IEC 61326-2-6; ТЕМ НЕ МЕНЕЕ ПЕРЕД НАЧАЛОМ ЭКСПЛУАТАЦИИ ПРИБОРА НЕОБХОДИМО ОЦЕНИТЬ ЭЛЕКТРОМАГНИТНЫЕ ХАРАКТЕРИСТИКИ ОКРУЖАЮЩЕЙ СРЕДЫ.

ДЛЯ ОБЕСПЕЧЕНИЯ НАДЛЕЖАЩЕГО ФУНКЦИОНИРОВАНИЯ ПРИБОРА ОПЕРАТОР ОБЯЗАН ПОДДЕРЖИВАТЬ СООТВЕТСТВИЕ ОКРУЖАЮЩЕЙ СРЕДЫ ТРЕБОВАНИЯМ ПО ЭЛЕКТРОМАГНИТНОЙ СОВМЕСТИМОСТИ.

ЗАПРЕЩАЕТСЯ ЭКСПЛУАТАЦИЯ ПРИБОРА ВБЛИЗИ ИСТОЧНИКОВ СИЛЬНЫХ ЭЛЕКТРОМАГНИТНЫХ ПОЛЕЙ (НАПРИМЕР, НЕЭКРАНИРОВАННЫХ ИСТОЧНИКОВ РАДИОЧАСТОТНОГО ИЗЛУЧЕНИЯ), КОТОРЫЕ МОГУТ НАРУШИТЬ РАБОТУ ПРИБОРА И ПРИВЕСТИ К НЕПРАВИЛЬНЫМ РЕЗУЛЬТАТАМ.





#### Общие сведения 2.

#### 2.1 Назначение

Фотометр INFINITE F50 PLUS — автоматизированный прибор с программным обеспечением Magellan, предназначенный для профессионального использования в лабораториях с целью измерения поглощения света (оптической плотности) в гомогенных жидких средах на 96-луночных микропланшетах в диагностике «in-vitro».

Он предназначен главным образом для выполнения диагностических анализов «in vitro» образцов тела человека, получаемых из выбираемого пользователем ферментного анализа (ELISA). Конкретная диагностическая информация и тип образца определяются выбранным анализом.

INFINITE F50 PLUS предназначен для измерения и оценки качественных, полуколичественных и количественных анализов в соответствии с запланированными диагностическими параметрами и техническими характеристиками прибора.

Изделие предназначено для профессионального использования в лабораториях обученным персоналом. Изделие не должно использоваться в домашних условиях или неквалифицированными лицами.





Примечание

В случае любой модификации фотометра INFINITE F50 PLUS или программного обеспечения Magellan, гарантия на прибор аннулируется, и он перестает соответствовать нормативным требованиям.

#### Примечание

При эксплуатации прибора для клинических диагностических исследований допускается использование только диагностических комплектов с маркировкой СЕ. Организация, эксплуатирующая прибор, должна удостовериться, что использование применяемого диагностического комплекта СЕ совместно с фотометром INFINITE F50 PLUS, имеющим маркировку СЕ, и его опциями, соответствует требованиям Директивы по диагностике in-vitro 2017/746 и другим относящимся государственным или местным предписаниям.

В случае использования фотометра INFINITE F50 PLUS в целях, отличных от представленных выше, или с программным обеспечением, отличным от Magellan, его соответствие требованиям IVD теряется, и ответственность за такое использование и необходимую аттестацию возлагается на организацию, эксплуатирующую прибор.



#### Примечание

Результаты, полученные с помощью фотометра INFINITE F50 PLUS, зависят от правильности эксплуатации прибора и микропланшетов, которая должна осуществляться в соответствии с инструкциями, приведенными в настоящем Руководстве, а также от используемых жидкостей (реагентов и других химических веществ). Необходимо строго соблюдать инструкции по работе и хранению прибора, а также инструкции по выполнению операций с образцами или реагентами. Следовательно, следует внимательно относиться к интерпретации полученных результатов.



Примечание Запрещается вскрывать корпус прибора. Данное действие влечет за собой потерю гарантии.

## 2.2 Принцип действия

Прибор INFINITE F50 PLUS представляет собой 8-канальный фотометр, предназначенный для измерения оптической плотности (OD) жидких сред на 96-луночных микропланшетах.

Для выбора длины волны при измерении абсорбции свет, создаваемый светодиодным источником, проходит через диск оптических фильтров.

Затем луч, разделенный восемью волоконными световодами, фокусируется на восемь лунок микропланшета. Прошедший через лунки свет направляется через восемь оптических линз на восемь фотодиодных детекторов.

Это означает, что INFINITE F50 PLUS способен одновременно измерить восемь лунок, т. е. один столбец 96-луночного микропланшета.

Для контроля и регулирования интенсивности излучения светодиодной лампы используется дополнительный волоконный световод (опорный канал), по которому свет проходит в обход микропланшета.

Полученные значения коэффициента пропускания преобразуются в значения оптической плотности (OD) согласно следующим формулам:

Коэффициент пропускания  $T = \frac{I}{I_o}$ 

I<sub>0</sub> = интенсивность падающего света

I = интенсивность детектированного света (после образца)

Оптическая плотность равна логарифму величины, обратной коэффициенту пропускания.

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



## 2.3 Профиль пользователя

# 2.3.1 Профессиональный пользователь - уровень администратора

Администратор должен иметь необходимое техническое образование, навыки и опыт, а также уметь распознавать и предотвращать опасности в процессе целевого использования изделия.

Администратор должен обладать обширными знаниями и уметь консультировать конечного или рядового пользователя по вопросам целевого использования изделия Тесап в задачах анализа.

Требуется знание прикладного программного обеспечения и хорошее владение английским языком.

## 2.3.2 Конечный или рядовой пользователь

Конечный или рядовой пользователь должен иметь необходимое техническое образование, навыки и опыт, а также уметь распознавать и предотвращать опасности в процессе целевого использования изделия.

Требуется знание прикладного программного обеспечения и хорошее владение национальным языком для данного объекта или английским языком.

## 2.3.3 Специалисты по ремонту

Специалист по ремонту должен иметь необходимое техническое образование, навыки и опыт, а также уметь распознавать и предотвращать опасности, связанные с ремонтом или обслуживанием изделия.

Требуется знание прикладного программного обеспечения и хорошее владение английским языком.



Примечание

Сведения о датах, продолжительности и частоте курсов обучения можно получить у вашего представителя службы поддержки покупателей.

Адрес и номер телефона приведены на веб-сайте <u>http://www.tecan.com/customersupport</u>



## 3. Приступая к работе

## 3.1 Распаковка и осмотр

## 3.1.1 Осмотр содержимого поставки

Комплект поставки:

- Внешний источник питания
- Шнур питания
- USB-кабель для подсоединения к внешнему компьютеру
- Руководство по использованию (PDF-файлы) на носителе данных
- USB-накопитель
  - о Программное обеспечение Magellan
  - Руководство по эксплуатации (PDF-файлы)
  - о Программные инструменты (например, Adobe Reader)





#### ПРЕДОСТЕРЕЖЕНИЕ

Примечание Во избежание потери данных и для предотвращения иных негативных последствий в случае вирусной атаки, запрещается снимать защиту записи с USB-накопителя.

ИСПЫТАНИЯ ПРИБОРА ПРОВОДИЛИСЬ С ИСПОЛЬЗОВАНИЕМ ПОСТАВЛЕННОГО USB-КАБЕЛЯ. ПРИ ИСПОЛЬЗОВАНИИ ДРУГОГО USB-КАБЕЛЯ ВОЗМОЖНО НЕШТАТНОЕ ФУНКЦИОНИРОВАНИЕ ПРИБОРА.

## 3.1.2 Порядок распаковки

- 1. Перед открыванием поставленной коробки осмотрите ее на предмет повреждений. Немедленно сообщите о любых обнаруженных повреждениях.
- Выберите место для установки прибора. Оно должно быть плоским и горизонтальным, не подверженным вибрации, защищенным от прямого солнечного света, свободным от пыли, паров растворителей и кислот. Расстояние между прибором и стеной должно быть не менее 5 см.
- 3. Извлеките прибор из коробки и поставьте его на выбранное место. Поднимать прибор следует осторожно.
- 4. Осмотрите прибор на предмет отсутствия плохо закрепленных, согнутых или сломанных деталей. Немедленно сообщите о любых обнаруженных повреждениях.
- 5. Сравните серийный номер прибора, указанный на его задней панели, с серийным номером в упаковочном листе. Немедленно сообщите в случае их расхождения.
- 6. Проверьте наличие принадлежностей по накладной.
- 7. Сохраните все упаковочные материалы, так как упаковка может понадобиться для транспортировки прибора в будущем.



## 3.2 Требования к питанию

Прибор автоматически определяет напряжение в сети питания, поэтому устанавливать нужное напряжение не требуется. Убедитесь, что напряжение питания имеет следующие характеристики:

Питание:	
Базовый прибор с адаптером пер. тока:	100–240 В, 50/60 Гц
Базовый прибор без адаптера пер. тока:	24 В пост. тока

Если данное напряжение не используется в вашей стране, обратитесь в местный сервисный центр Tecan.

Прибор допускается подключать только к электросети с защитным заземлением.



ПРЕДОСТЕРЕЖЕНИЕ ЗАПРЕЩАЕТСЯ ПОДСОЕДИНЯТЬ ПРИБОР К ЭЛЕКТРОСЕТИ С ДРУГИМ ДИАПАЗОНОМ НАПРЯЖЕНИЙ. ВКЛЮЧЕНИЕ ПРИБОРА, ПОДСОЕДИНЕННОГО К ТАКОЙ ЭЛЕКТРОСЕТИ, ПРИВЕДЕТ К ЕГО ПОВРЕЖДЕНИЮ.



## ПРЕДОСТЕРЕЖЕНИЕ

НЕ ИСПОЛЬЗУЙТЕ ВМЕСТО ОТДЕЛЯЕМЫХ ШНУРОВ ПИТАНИЯ ДРУГИЕ ШНУРЫ С НЕСООТВЕТСТВУЮЩИМИ ХАРАКТЕРИСТИКАМИ.

## 3.3 Требования к окружающей среде

Прибор следует устанавливать на плоскую горизонтальную поверхность, свободную от пыли, паров растворителей и кислот.

Для обеспечения точности полученных результатов необходимо исключить воздействие вибрации и прямого солнечного света.

Температура окружающего воздуха:	
Работа	от 15 до 35 °C (от 59 до 95 °F)
Хранение	от -30 до 60 °С (от -22 до 140 °F)
Относительная влажность:	от 20 до 80 %, отсутствие конденсации при рабочей температуре



# 3.4 Требования к системе

	Минимальные	Рекомендуемые
пк	Windows-совместимый ПК с процессором типа Pentium или совместимым с ним, работающим на частоте 1 ГГц	2 ГГц (Dual Core)
Операционная	Windows 10 (32-разрядная)	
система	Windows 10 (64-разрядная) Версия: Pro	
Память	Windows 10 (32-разрядная): 1 Гб (ОЗУ)	2 Гб (ОЗУ)
	Windows 10 (64-разрядная): 2 Гб (ОЗУ)	4 Гб (ОЗУ)
Свободное место на жестком диске	3 Гб	5 Гб
Монитор	Графика Super VGA	
Разрешение	1024 х 600 и выше	1920 x 1080
Число цветов	256	
Мышь	Мышь Microsoft или совместимое указывающее устройство	
Интерфейс	1 x USB 2.0	2 x USB 2.0 1 x RS232 (последовательный)
Устройства	Windows 10:	
	Видеокарта с поддержкой DirectX 9 и драйвером WDDM 1.0 или более поздней версии	
.NET	Microsoft .NET Framework 3.5: В Windows 10, если .NET Framework (3.5) еще не установлена, поступит запрос на ее установку.	
Windows Installer	<ol> <li>3.1</li> <li>Если эта версия не установлена, программа установки/обновления установит ее.</li> </ol>	
Microsoft Excel	2007, 2010, 2013, 2016 (32-разрядная), 2019 (32-разрядная), Microsoft Excel 365 (32-разрядная) Поддерживаются только 32-разрядные версии! Начальные версии НЕ поддерживаются!	2010 (32-разрядная), 2019 (32-разрядная)



## 3.5 Включение прибора

В этом разделе приведены действия, которые необходимо выполнить перед включением прибора.



#### ПРЕДОСТЕРЕЖЕНИЕ ДЛЯ УСТРАНЕНИЯ ОПАСНОСТИ ВЫЗВАННОГО КОНДЕНСАЦИЕЙ КОРОТКОГО ЗАМЫКАНИЯ, ПЕРЕД УСТАНОВКОЙ И ВКЛЮЧЕНИЕМ ПРИБОРА ЕГО НЕОБХОДИМО ВЫДЕРЖАТЬ В ТРЕБУЕМЫХ УСЛОВИЯХ ОКРУЖАЮЩЕЙ СРЕДЫ В ТЕЧЕНИЕ НЕ МЕНЕЕ ТРЕХ ЧАСОВ.

После выполнения приведенных выше требований можно приступать к установке, выполняемой следующим образом:

- 1. Подсоедините прибор к внешнему компьютеру через USB-кабель.
- 2. Убедитесь, что выключатель питания на левой панели находится в выключенном состоянии.
- 3. Установите шнур питания в гнездо питания на левой панели.
- 4. Включите прибор выключателем питания на левой панели.

Прибор будет готов к измерению микропланшетов после установки программного обеспечения.

#### Примечание

Перед началом измерений необходимо убедиться, что позиция А1 микропланшета находится в надлежащем положении.

Измерения микропланшетов возможны только без крышек.

Во избежание искажения результатов окружающим светом перед началом измерения закройте крышку транспорта планшета.



Примечание Всегда извлекайте микропланшет из фотометра сразу после завершения измерения.





#### Программное обеспечение 3.6

#### 3.6.1 Введение / Назначение

Программное обеспечение Magellan, используемых для управления фотометром и обработки данных, входит в комплект поставки прибора.

Magellan представляет собой универсальное программное обеспечение для управления прибором и обработки данных, получаемых при проведении измерений на микропланшетах с использованием измерительного оборудования компании Tecan.

Поставляются две различные версии программного обеспечения Magellan:

- Magellan и
- **Magellan Tracker**

Magellan — версия программного обеспечения, предназначенная для измерений конечной точки, кинетических анализов и анализов с несколькими метками на приборе INFINITE F50 PLUS в соответствии с назначением; см. раздел 2.1 Назначение.

Magellan Tracker — версия программного обеспечения, обеспечивающая, в дополнение к функциональным возможностям Magellan, соблюдение требований части 11 положения 21 Свода федеральных нормативных актов США, разработанного Управлением по контролю за продуктами и лекарствами США (FDA 21 CFR part 11).

# Примечание

Важно понимать, что правильная установка оборудования и программного обеспечения Magellan сама по себе не обеспечивает соответствия всем применимым законам и требованиям. Необходимо также установить соответствующие правила в отношении процессов и стандартных рабочих процедур, включая подтверждение достоверности и контроль качества.

#### 3.6.2 Порядок установки

Для установки программного обеспечения вставьте USB-накопитель в USBпорт и выполните следующие действия:

- 1. Автоматически запустится мастер установки Magellan, который поможет выполнить процедуру установки. Если мастер не запустился автоматически, запустите программу E:\Tecan.exe (где «Е» — буква диска USB-накопителя).
- 2. Для запуска процедуры установки выберите 'Magellan software' и (в зависимости от приобретенной версии) 'Install Magellan' или 'Install Magellan Tracker', затем следуйте подсказкам мастера.
- 3. Нажмите кнопку Установить для запуска процедуры установки программного обеспечения.
- 4. Для продолжения щелкните в поле I accept the terms of the license agreement и нажмите кнопку Next.
- 5. На открывшейся странице Сведения о клиенте введите имя пользователя и название организации.
- 6. На странице Конфигурация выберите язык.



#### 3. Приступая к работе



- 7. Страница Use for regulated environments: для продолжения нажмите кнопку Next
- 8. Нажмите кнопку Установить для запуска процедуры установки.
- 9. Нажмите кнопку Готово для завершения установки и выхода из программы установки.

Программное обеспечение можно запустить через меню **Пуск** Windows путем выбора **Magellan** в группе программ **Tecan**.





Примечание Совершенно необходимо, чтобы пользователь, выполняющий установку программного обеспечения, имел права администратора на данном компьютере.

Примечание Magellan V7.5 невозможно установить параллельно с другими версиями Magellan.

По умолчанию, все файлы, относящиеся к программному обеспечению Magellan for Infinite F50, хранятся в подкаталогах каталога:

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

## 3.6.3 Проверка установки

Выполните проверку правильности установки программного обеспечения Magellan при помощи программы автоматической проверки установки:

Запустите программу *TecanIQ.exe* из каталога установки по умолчанию (C:\*Program Files*\*TecanWagellan*) или через меню **Пуск: Start > Programs > Tecan > MagellanIQ**.

Нажмите кнопку **Проверить**, чтобы начать проверку установки. Все установленные компоненты должны иметь статус **ОК**. При наличии сообщения о какой-либо потенциальной проблеме обратитесь к местному торговому представителю Tecan.

Для закрытия программы проверки установки нажмите кнопку **Отмена** или **Выход**.



Примечание

Процедуру проверки установки следует повторять каждый раз после установки программного обеспечения Magellan или обновления до более новой версии.



#### 3.6.4 Начало работы с программным обеспечением Magellan

Основным видом пользовательского интерфейса в программном обеспечении Magellan являются мастера.

Стандартные мастера в Magellan представляют собой рабочие модули с пошаговыми инструкциями для выполнения сложных процедур.

Для некоторых операций имеются меню в головной панели. **Меню** обеспечивает традиционный способ использования программного обеспечения: выбор требуемого пункта меню из главного меню. Все последующие действия запускаются немедленно, или же отображается диалоговое окно, позволяющее сделать дальнейший выбор или ввести данные.

#### Пользовательский интерфейс – Список мастеров

После загрузки программного обеспечения Magellan отображается Список мастеров.

Список мастеров - Magellan Tracker	
$\frown$	Выберите одну из следующих
Запуск измерения	Мастер "Начало измерения" поможет Вам провести измерение.
Оценка результатов	<ul> <li>Вы можете использовать метод или получить необработанные данные</li> </ul>
Подписать	•
Созд./редактир. списка идент. образцов	•
Созд./редактир. метода	•
/	
Выход Magellan	

Чтобы запустить какой-либо мастер, дважды щелкните по нему или

выберите его и нажмите кнопку Далее

#### Мастер Запуск измерения

В мастере Запуск измерения предусмотрены следующие параметры:

- Получить исходные данные используется для быстрого и простого получения исходных данных путем установки требуемых параметров измерения и запуска измерения.
- Использование предопределенного метода применяется для выполнения измерений согласно ранее определенным методам.
- Запуск избранного метода используется для выбора одного из наиболее часто используемых методов из списка нумерованных значков.

По окончании измерения будет создан файл рабочего пространства.



#### Мастер Оценка результатов

Мастер **Оценка результатов** используется для просмотра исходных данных и оценки результатов. Также возможен просмотр параметров оценки данных и повторная оценка данных.

#### Мастер Прикрепление подписи

Мастер **Прикрепление подписи** используется для подписи файлов методов и рабочих пространств. Данная функция доступна только в Magellan Tracker.

#### Мастер Создание/редактирование списка идентификаторов образцов

Мастер **Создание/редактирование списка идентификаторов образцов** используется для создания новых и редактирования существующих списков идентификаторов образцов.

#### Мастер Создать/редактировать метод

Мастер Создание / редактирование метода используется для определения и редактирования методов.



Примечание Более подробную информацию по использованию Magellan см. в «Руководстве по использованию» программного обеспечения Magellan.

Следует иметь в виду, что некоторые функции, описанные в «Руководстве по использованию» программного обеспечения Magellan, могут оказаться недействительными в комбинации Magellan V7.5 с INFINITE F50 Plus. Тем не менее, в этом «Руководстве по использованию» представлены все необходимые сведения.



Примечание Пример анализа ELISA подробно рассматривается в главе 6 Пример применения.

Файлы метода Sunrise, созданные при помощи программного обеспечения Magellan V7.x или более нового, открываются с помощью Magellan V7.5, при этом параметры измерения конвертируются автоматически.

Методы Sunrise, созданные в более ранних версиях, необходимо конвертировать. Для этого в меню «Дополнительно/Обработка файла» выберите пункт «Конвертировать из».



## 3.7 Magellan – Редактор параметров измерения

Редактор параметров измерения используется для настройки рабочих процессов. Для создания рабочего процесса достаточно перетащить в последовательность шаги, выбранные в соответствии с приложением. Затем рабочий процесс приложения можно просмотреть на панели рабочего процесса. Каждый шаг рабочего процесса (программный элемент) можно скопировать и вставить (используя стандартные комбинации клавиш Windows Ctrl+C и Ctrl+V или контекстное меню мыши) или переместить в выбранное положение.

Тараметры измерения - infinite F50			×
🔶 Лаб. оборудование 🔹	🔷 🔻 Планшет	1 Выбор	
🔶 Планшет 🍝 Часть планшета	Onpeg. nnawueta: [GRE96R] - Greiner 96 Flat Transparent	Детали Использ. часть планшета	ine bogesterio
У Изнерения 🛞 Фолтическая плотность	№ № Часть планиета А1:H12	2	
Действия 🔅 Встряхивание Перенести планшет	Aness pone Weepervee: 445 mm V Transe: 445 mm V Vee: Mencal V		
<ul> <li>Кинетическое измерение</li> <li>Кинетической цикл</li> <li>Условия кинетики</li> </ul>			
Дополнительно 🙁 Кончентарии 🔶 Загрос пользователя Э Ожидать (тайкер) Э Инкубация			
Панель управления	Панель рабочего процесса	Информа панель	ационн
			×
Число планшетов: 1 🔹 Справка Отнена <<<Назад		Выберите параметры измерения	

Редактор параметров измерения состоит из следующих элементов, подробно описываемых далее:

- Панель управления
- Панель рабочего процесса
- Информационная панель



## 3.7.1 Панель управления

**Панель управления** разделена на пять частей. Каждая часть содержит программные элементы, используемые для создания различных рабочих процессов.

Для создания рабочего процесса необходимо дважды щелкнуть выбранный программный элемент или перетащить его на панель рабочего процесса.

При использовании фотометра INFINITE F50 PLUS доступны следующие программные элементы:

Лабораторное оборудование	Планшет Часть планшета
Измерения	Оптическая плотность
Действия	Встряхивание Перенести планшет
Кинетика	Кинетический цикл Условия кинетики
Дополнительно	Комментарии Запрос пользователя Ожидать (Время) Инкубация

#### Лабораторное оборудование

#### Планшет

Программный элемент **Планшет** используется для выбора формата планшета из раскрывающегося списка **Определение планшета**. Для просмотра дополнительной информации по выбранному планшету нажмите кнопку **Детали**.

🔷 🔻 Планшет		1
Опред. планшета:	[CO596ft] - Corning 96 Flat Transparent	🖌 Детали
		Использ, часть планшета

#### Часть планшета

По умолчанию программный элемент **Часть планшета** находится в свернутом состоянии. В раскрытом состоянии на нем отображается 96луночный микропланшет. Для измерения отдельных лунок щелкните по выбранной лунке. Для измерения диапазона лунок обведите выбранные лунки рамкой путем перетаскивания ее границ. Для предварительного просмотра планшета в увеличенном масштабе нажмите кнопку **Детали**.

**		
1       2       3       4       5       6       7       8       9       10       11       12         B       Image: Constraint of the state	Детали	



#### Независимые части планшета

Предусмотрена возможность выбора независимых частей планшета:



Чтобы указать второй диапазон лунок, нажмите клавишу **Ctrl** и обведите рамкой выбранные лунки.

#### Измерения

#### Оптическая плотность

Программный элемент **Оптическая плотность** используется для выполнения измерений оптической плотности. Введите или выберите соответствующие параметры.

В двух раскрывающихся списках отображаются, в соответствии с установленными фильтрами поглощения, доступные длины волн измерения и референсные длины волн. Если фильтры не определены, раскрывающиеся списки будут пусты.

🍦 🔻 Оптич	еская плотнос	гь		3
Длина волны Измерение:	492 nm	~	Метка Имя: Метка1	
🔲 Эталон:	405 nm	~		

#### Действия

#### Встряхивание

Программный элемент Встряхивание используется, когда перед началом измерения или между кинетическими циклами планшет необходимо встряхнуть.

📉 🔻 Встрякивание	4
Параметр Продолжит: 1 📚 с Интенсивность: Wide 💌 Амплитуда: 14,1 mm Частота: 2,1 Hz	W ALTE DOW OF YMA

Введите требуемые параметры:

Продолжит.	Введите продолжительность процесса встряхивания.
Интенсивность	Введите требуемый режим встряхивания. При выборе режима встряхивания отображаются амплитуда и частота выбранного режима.



#### 3. Приступая к работе

Предусмотренные режимы встряхивания приведены в главе 4.1.1 Встряхивание микропланшета.

Чтобы вставить новый программный элемент, нажмите ссылку <u>Ждите пару</u> секунд. Подробное описание см. на стр. 25.

#### Кинетика

#### Кинетический цикл

Программный элемент Кинетический цикл используется для выполнения нескольких последовательных измерений, которые могут осуществляться через заданные интервалы времени.

😡 🔻 Кинетический цикл		3
Циклы () Кол-во циклов: 2 () Продолжит.	Кинетический интервал ✓ Используйте кинетический интервал: ④ Время: 00:01:00 ♀ (чч:ми:сс) ○ Время: 60000 ♀ мс	

Введите требуемые параметры:

Циклы	Количество циклов: введите число или выберите с помощью кнопок со стрелками вверх или вниз количество шагов фактических измерений (2 – 1000). Продолжит.: введите продолжительность цикла в формате чч:мм:сс.
Кинетический интервал	Используйте кинетический интервал: введите интервал времени (чч:мм:сс или мс).

#### Условия кинетики

Программный элемент **Условия кинетики** используется для определения действий, выполняемых в конкретном цикле.



Если ввести значение **2** в поле **Выполнить команды при цикле** процесса кинетического измерения, в котором содержится, например, шаг **Встряхивание**, встряхивание будет выполняться только на шаге 2.

#### Примечание

Кинетические условия, такие как встряхивание, следует вставить сразу после программного элемента «Кинетический цикл», с тем чтобы обеспечить оптимальный результат воспроизводимости.

Для получения сопоставимых результатов пользователям рекомендуется создать соответствующие сценарии до проведения измерений и использовать один и тот же сценарий для всех аналогичных кинетических измерений.





#### Дополнительно

#### Комментарии

Программный элемент Комментарии используется для ввода в текстовом поле замечания или пояснения к текущему измерению.

🚛 🔻 Комментарии	6
Комментарии:	

#### Запрос пользователя

Программный элемент **Запрос пользователя** сообщает оператору о необходимости выполнения в определенное время рабочего процесса какого-либо действия.

🚯 🔻 Запрос пол	ьзователя	7
Текст:		

Если, например, для извлечения планшета используется программный элемент **Перенести планшет**, то введенный текст сообщает оператору о необходимости выполнения требуемых действий. Это сообщение отображается в диалоговом окне, где для продолжения процесса измерения следует нажать кнопку **OK**.

#### Ожидать (Время)

Программный элемент **Ожидать (Время)** используется для задания определенного времени ожидания до начала выполнения следующего шага процесса.

#### Введите требуемое время в поле Время ожидания.

🧐 🔻 Ожидать (таймер)		5
Таймер Время ожидания: 00:01:00 🗘 (чч:мм:сс)	Параметры Ожидание для впрыска Игнорир. ожидание посл. кинет. цикла	

Введите требуемые параметры:

Таймер	Введите Время ожидания (чч: мм: сс)	
Параметры	Игнорир. ожидание посл. кинет. цикла: Если программн шаг Ожидать (Время) является последним действием в ходе кинетического анализа, в последнем цикле он буде игнорироваться.	

#### Инкубация

🧐 🔻 Инкубация	4
Таймер Время инкубации: 00:01:00 文 (чч:мм:сс)	

Введите требуемые параметры инкубации:

Время	Введите общее время (не менее 5 с)
инкубации	



## 3.7.2 Панель рабочего процесса

Панель рабочего процесса представляет собой окно, в котором отображается сценарий измерения и можно определить и изменить его параметры.

Существует два способа вставки программного элемента из Панели управления в Панель рабочего процесса:

- Выберите программный элемент на Панели управления двойным щелчком. Выбранный элемент вставляется в Панель рабочего процесса сразу после предыдущего программного элемента.
- Выберите программный элемент на **Панели управления** и перетащите его на **Панель рабочего процесса** в требуемое положение.

Нумерация программных элементов осуществляется в соответствии с их местом в последовательности.

После помещения программного элемента на Панель рабочего процесса его настройки и параметры становятся доступными для изменения.

Отдельные программные элементы на **Панели рабочего процесса** можно свернуть для отображения более важной информации или развернуть для получения доступа ко всем изменяемым функциям. Для переключения между двумя режимами просмотра нажмите один из треугольников или , расположенных рядом с названием программного элемента.

По умолчанию, сначала на **Панели рабочего процесса** в редакторе параметров измерения отображается программный элемент **Планшет**, затем элементы **Часть планшета** (в свернутом виде) и **Оптическая плотность**.

Текущие выбранные программные элементы обозначаются на **Панели рабочего процесса** желтой линией на верхней границе.

Если программный элемент содержит ошибки или недействителен в текущем процессе, он помечается флагом ошибки, а его номер отображается красным цветом. Подробные сведения об ошибки отображаются на **Информационной панели**. Если в рабочем процессе имеются ошибки, выбор параметров измерения невозможен.

#### Иерархия элементов

В Панели рабочего процесса действует следующая иерархия элементов:

- 1. Планшет
- 2. Часть планшета (Диапазон)

Любой шаг измерения можно вставить сразу после элементов «Планшет» или «Диапазон». Изменить последовательность выполнения отдельного стрипового компонента можно с помощью функций **Разблокировать** и **Отступ**. Укажите элемент на **Панели рабочего процесса**, нажмите правую кнопку мыши и выберите пункт **Разблокировать** или **Отступ**.

Другие элементы **Панели управления** можно вставить в иерархию рабочего процесса следующим образом:

Первый элемент **Диапазон** вставляется сразу после элемента **Планшет**; затем вставляются все последующие элементы **Диапазон**.

Кинетические шаги можно задавать в элементах Планшет или Диапазон.

Шаги Запрос пользователя, Комментарии и Ожидать можно задавать в элементах Планшет или Диапазон.



#### 3.7.3 Информационная панель

На расположенной в правой части экрана **Информационной панели** отображаются сведения по выбранному в данный момент программному элементу. Здесь также отображаются все предупреждения и ошибки.

## 3.8 Magellan - Определение измерений

В этой главе представлены некоторые примеры определения параметров различных измерений.

#### 3.8.1 Определение измерений конечной точки

В этом примере описывается Измерение конечной точки для оптической плотности во всех лунках 96-луночного микропланшета.

- 1. Выберите 96-луночный микропланшет в раскрывающемся списке Описание планшета.
- 2. По умолчанию, для измерения выбраны все лунки 96-луночного микропланшета.
- 3. Введите требуемые длины волн измерения и референсные длины волн.

🔷 🔻 Планшет			1
Опред. планшета: [О	OS96ft] - Corning 96 Flat Transparent		<ul> <li>Детали</li> <li>Использ. часть планшета</li> </ul>
<ul> <li>Часть п</li> <li>1 2 3 4</li> <li>В</li> <li>В</li></ul>	ланиета 5 6 7 8 9 10 11 12 6 6 7 8 9 0 10 11 12 6 6 7 8 9 10 10 10 10 7 7 7 7 7 7 7 7 7 7 7 7 7 7 7 7 7 7 7	אתנ	2
С О О О О О О О О О О О О О О О О О О О	IT/IY4ECKA3 ПЛЮТНОСТЬ ПНЫ е: 450 nm ▼ 4: 520 nm ▼	Метка Имя: Метка1 💌	3



## 3.8.2 Определение измерений с несколькими метками

В этом примере описывается **Измерение с несколькими метками для** оптической плотности в заданном диапазоне (A1:E7) 96-луночного микропланшета. Предполагается измерить три метки оптической плотности.

- 1. Выберите 96-луночный микропланшет в раскрывающемся списке Описание планшета.
- 2. По умолчанию, для измерения выбраны все лунки 96-луночного микропланшета.

Нажмите **•**, чтобы развернуть элемент **Часть планшета**. Затем выберите требуемый диапазон (A1:E7).

- 3. Введите требуемую длину волны измерения.
- 4. Вставьте два дополнительных элемента Оптическая плотность и введите для них длины волн измерения.

Chype, Innakusers       LC0598tj - Coming 96 Flat Transparent       L       LC0598tj - Coming 96 Flat Transparent         Image: Comparison of the comparison of t					
Vacrь планиета Vacrь планиет	Опред. планшета:	[COS96ft] - Corning 96 Flat Transparent			🖌 Детали
Vactor Intelliere   I					Использ, часть план.
A 5 6 7 8 9 10 11 12       Arram         B 6       A 5 6 7 8 9 10 11 12         C 7 00000000000000000000000000000000000	🏾 🔞 🔻 4ac	гь планшета			
Image: Antriveckas Informets     Anisha Bonkel     Image: Distribute Bonkel </th <th>A 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0</th> <th></th> <th>Детали</th> <th></th> <th></th>	A 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0		Детали		
Измерение:       405 nm         Эталон:       405 nm         Иня:       Метка         Иня:       Метка         Измерение:       492 nm         Эталон:       405 nm         Иня:       Метка         Иня:       Метка         Иня:       Метка         Измерение:       492 nm         Эталон:       405 nm         Мина:       Метка         Иня:       Метка         Иня:       Метка         Измерение:       620 nm         Эталон:       405 nm		Оптическая плотность			
Эталон:     405 nm       Линна Волны:     492 nm       Измерение:     492 nm       Эталон:     405 nm	Измер	а волны рение: 405 nm 🗸	Метка-	Метка1	
Изирение:       492 пл         Эталон:       405 пл         У       Оптическая плотность         Изиврение:       620 пл         Уня::       Метка         Изиврение:       620 пл         Эталон:       405 пл		алон: 405 nm 🗸			
Длина волны       Чог пл       Чика:       Мика:					
Изнерение: <u>492 nm</u> Эталон: <u>405 nm</u> Иня: <u>Метка2</u> Иня: <u>Метка2</u> Иня: <u>Метка3</u> Иня: <u>Метка3</u> Эталон: <u>405 nm</u>		Оптическая плотность			
Эталон:     405 nm       ↓     •       Длина волны       Изнерение:       620 nm       Эталон:       405 nm	4 -	з волны	Метка	]	
Дляна волны         Измерение:         620 пт           Эталон:         405 пт         Иня:	Длин- Измер	оптическая плотность а волны ение: 492 nm 💌	Метка- Имя:	Метка2	-
Дляна волны         Калина волны         Иня         Метка           Изиврение:         620 nm         Иня:         Метка         Иня:         Метка           Эталон:         405 nm         Иня:         Метка         Метка         Метка         Метка	Длин- Измер Эт	а волны ение: 492 nm v алон: 405 nm v	Иня:	Метка2	
Изнерение: <u>620 nm</u> Эталон: <u>405 nm</u> Иня: <u>Метка3</u>	Длин- Измер Эт	иптическая плотность а волны нение: 492 nm апон: 405 nm Оптическая плотность	Иня:	Метка2	
Эталон: 405 nm	Длин- Измер Эт Длин-	иптическая плотность а волны нение: 492 nm ♥ алон: 405 nm ♥ Оптическая плотность а волны	Иня:	Метка2	
	Алин. Измер Эт Алин. Измер	Оптическая плотность     аволны     аволны     492 nm     405 nm     Оптическая плотность     волны     волны     волны     векие: 620 nm     С	Иня:	Метка2 💌	
	Длин- Измер Эт Измер Эт	ОПТИЧЕСКАЯ ПЛОТНОСТЬ     а волны     405 nm     Ф     ОПТИЧЕСКАЯ ПЛОТНОСТЬ     В волны     волны     620 nm     Ф     Оптическая плотность	Метка Иня: Иня: Иня:	Метка2 v	



## 3.8.3 Определение кинетических измерений

В этом примере описывается кинетическое измерение 96-луночного микропланшета.

- 1. Выберите 96-луночный микропланшет в раскрывающемся списке Описание планшета.
- 2. Вставьте программный элемент **Кинетический цикл** между элементами «Часть планшета» и «Оптическая плотность».
- 3. Циклы/Количество циклов: 50
- 4. Кинетический интервал (интервал между измерениями): выберите **Используйте кинетический интервал** и введите: 2 минуты 30 секунд.
- 5. Задайте элемент Оптическая плотность путем ввода требуемой длины волны измерения.

🔶 🔻 Планшет	1
Опред. планшета:	[COS96ft] - Corning 96 Flat Transparent   Kcones, vacte nnarwera
🔹 🔻 Част	ъпланшета 2
1       2       3         B       •       •       •         C       •       •       •         D       •       •       •         F       •       •       •         G       •       •       •         H       •       •       •	A 5 6 7 8 9 10 11 12 Aerank
<b>•</b>	Кинетический цикл 3
Цикль () Кол Про	н л-во циклов: 50 ♀ одолжит. 50 ♀ Одолжит.
	Оптическая плотность
	Длина волны Измерение: 492 nm Эталон: 405 nm
	- 03



## 3.8.4 Создание отступов и разблокирование программных элементов

Создание отступов и разблокирование программных элементов позволяет изменить рабочий процесс в ходе измерений.

Операции, предусмотренные во всех программных элементах с одинаковым отступом, выполняются последовательно. Единственной зависимостью между этими элементами является то, что следующее действие начинается сразу после завершения предыдущего.

Больший отступ одного элемента относительно другого указывает на зависимость между ними. Это означает, что действие параметров, определенных в первом элементе, распространяется также и на второй (с отступом) элемент.

В следующем примере представлен процесс определения **Меток** кинетических измерений для двух **Меток оптической плотности**. На этом примере видно, что два элемента **Оптическая плотность** зависят от элемента **Кинетический цикл**, который, в свою очередь, зависит от элемента **Часть планшета**, зависящего от элемента **Планшет**. Задайте следующие параметры для этого примера:

- 1. Планшет: например, Грейнер, 96, плоский, прозрачный
- 2. Кинетический цикл/Количество циклов: 5
- 3. Оптическая плотность/длина волны, метка 1: 450 нм
- 4. Оптическая плотность/длина волны, метка 2: 492 нм

Теперь Панель рабочего процесса отображается в следующем виде:

🔷 🔻 Планшет	1
Опред. планшета: [GRE96ft] - Greiner 96 Flat Transparent	Детали Использ. часть планшета
🎪 🔻 Часть планшета	2
1       2       3       4       5       6       7       8       9       10       11       12         B       Image: Constraint of the state	
	3
Циклы © Кол-во циклов: 5 Продолжит.	
👌 🔻 Оптическая плотность	4
Длина волны Измерение: 450 nm ✓ □ Эталон: 405 nm ✓	
🔻 Оптическая плотность	5
Алина волны Измерение: 492 nm Эталон: 405 nm Умя: Метка2	
	203

Такое определение параметров задает следующий рабочий процесс:



сначала измеряется поглощение во всех лунках 96-луночного микропланшета на длине волны 450 нм, затем на длине волны 492 нм. Оба измерения оптической плотности выполняются в ходе пяти кинетических циклов.

Создание отступа второго элемента Оптическая плотность до одного уровня с элементом Кинетический цикл приводит к изменению рабочего процесса. Выберите второй элемент Оптическая плотность и нажмите правую кнопку мыши. Выберите в контекстном меню пункт Разблокировать стрип. Теперь Панель рабочего процесса отображается в следующем виде:

_
ר
ета
:
;
2

Согласно этому рабочему процессу сначала выполняется **Кинетическое** измерение оптической плотности за пять циклов на длине волны 450 нм. По его завершении выполняется **Измерение конечной точки оптической** плотности на длине волны 492 нм.

## 3.9 Оптимизация эксплуатационных характеристик

Прибор прошел в полном объеме заводские испытания на соответствие заявленным характеристикам (подробнее см. 4.4.2 Характеристики измерения).

Максимальная точность измерений достигается при соблюдении приведенных ниже рекомендаций.

## 3.9.1 Расположение прибора

Место установки прибора должно отвечать требованиям, представленным в разделе 3.3 Требования к окружающей среде.



## 3.9.2 Процедура эксплуатации

#### Общие сведения

- Рекомендуется придерживаться стандартных рабочих процедур анализа.
- Наилучшая воспроизводимость достигается при использовании длины волны измерения, равной длине волны максимальной оптической плотности данного раствора.
   Важно использовать длину волны максимальной оптической плотности, если кривая поглощения образца лежит в узком диапазоне длин волн.

Имейте в виду, что проведение измерений на наклонной части пика оптической плотности ограничивает точность полученных значений.

- После измерения каждого микропланшета, ознакомьтесь с информацией по процедуре проверки, которая приводится на упаковке диагностического комплекта.
- Используйте фильтры поглощения, рекомендованные для фотометра INFINITE F50 PLUS.

#### Микропланшеты

- Допустимые типы микропланшетов для данного прибора указаны в главе 4.4.3 Микропланшеты.
- Наилучшие результаты достигаются при использовании микропланшетов с плоским дном. Тип микропланшета может влиять на результаты измерения.
   Особую осторожность соблюдайте при использовании планшетов с лунками с С-, U- или V-образным дном или лунками в стрипах, так как результаты измерений могут несколько отличаться от описанных в спецификациях, приведенных в этом документе. Тип микропланшета, используемого с фотометром INFINITE F50 PLUS, должен соответствовать применению.
- Используйте только совершенно чистые микропланшеты.
- Не допускайте оседания пыли на растворах или на микропланшете, если микропланшет оставляют стоять в течение некоторого времени перед измерением.
   Каждый раз, когда микропланшет находится вне прибора, его рекомендуется закрывать защитной крышкой.
- Неточности в количестве раствора в лунках сильно влияют на полученные результаты, если используются малые количества раствора.
- Форма мениска раствора может вызвать неточности в результатах, особенно при использовании малых количеств раствора



# 4. Описание прибора

## 4.1 Описание прибора

В фотометре INFINITE F50 PLUS предусмотрены следующие режимы измерения оптической плотности:

измерение конечной точки, кинетическое измерение и измерение с несколькими метками.

### 4.1.1 Встряхивание микропланшета

Фотометр INFINITE F50 PLUS позволяет встряхивать микропланшет перед началом измерений. Встряхивание микропланшета также может выполняться между циклами кинетического измерения.

Настройка режима встряхивания выполняется в Magellan.



## ПРЕДОСТЕРЕЖЕНИЕ

ПРИМЕНЕНИЕ РЕЖИМА ВСТРЯХИВАНИЯ ПРИ ВЫСОКОМ УРОВНЕ ЖИДКОСТИ В ЛУНКАХ МОЖЕТ ПРИВЕСТИ К ЕЕ ПРОЛИВАНИЮ.

Режимы встряхивания INFINITE F50 PLUS:

Режим встряхивания	Амплитуда встряхивания	Частота встряхивания
ИНТЕНСИВНЫЙ	2,8 мм	14,3 Гц
ОБЫЧНЫЙ	4,4 мм	9,2 Гц
СЛАБЫЙ	4,4 мм	7,8 Гц
ШИРОКИЙ	14,1 мм	2,0 Гц

## 4.2 Описание прибора

На рисунке ниже показаны компоненты прибора.



2023-06



#### 4. Описание прибора

Индикатор состояния прибора имеет следующие режимы:

- Мигающий зеленый: прибор не подключен к Magellan;
- Зеленый: прибор подключен и готов к выполнению измерения;
- Красный: выполняется измерение.

На левой панели прибора размещены USB-порт, выключатель и гнездо питания.



Паспортная табличка прикреплена к днищу прибора.

Образец паспортной таблички



Содержание паспортной таблички (например, название модели и номер изделия) может зависеть от конкретной модели.

Перечень различных приборов, на которые распространяется настоящее Руководство по эксплуатации, приведен в Заявлении о соответствии ("Declaration of Conformity") на последней странице данного документа.



## 4.3 Описание диска фильтров

В стандартном диске фильтров, поставляемом в комплекте фотометра INFINITE F50 PLUS, установлены четыре узкополосных интерференционных фильтра с фиксированной длиной волны (405, 450, 620 и 492 нм). Всего в диск можно установить до восьми фильтров. Для заказа дополнительных фильтров обратитесь к местному торговому представителю Tecan.

В стандартном диске фильтры установлены в следующем порядке:

Позиция фильтра	Длина волны фильтра
1	405 нм
2	450 нм
3	620 нм
4	492 нм
5 - 8	Пустые позиции для фильтров



После выбора длины волны измерения диск передвигается таким образом, чтобы луч света проходил через требуемый фильтр.



Примечание Подробнее об определении нового фильтра см. раздел 7.5.2 Определение фильтров.

## 4.4 Технические характеристики прибора

В следующей таблице приведены технические характеристики фотометра INFINITE F50 PLUS.

## 4.4.1 Общие характеристики

ΠΑΡΑΜΕΤΡ	ЗНАЧЕНИЕ	
Питание	Питание:	
Внешний источник питания	Базовый прибор с адаптером пер. тока: 100–240 В, 50/60 Гц, макс. 1,2 А (автоопределение, категория перенапряжения II)	
	Базовый прибор без адаптера пер. тока: 24 В пост. тока (категория перенапряжения I)	
Потребляемая мощность INFINITE F50 PLUS	Режим ожидания: прибл. 12 ВА Рабочий режим: макс. 30 ВА	
Габаритные размеры	Ширина: 34,7 см (13,66 дюйма) Глубина: 18,9 см (7,44 дюйма) Высота: 13,4 см (5,28 дюйма)	
Масса	2,6 кг (включая источник питания)	
Температура окружающе	его воздуха:	
Работа	от 15 до 35 °C (от 59 до 95 °F)	
Хранение	от -30 до 60 °C (от -22 до 140 °F)	
Относительная влажность	от 20 до 80 %	
Степень загрязнения	2	
Метод утилизации	Загрязненные отходы	
Экологические требования	Более подробную информацию см. в разделе 3.3 Требования к окружающей среде.	


## 4.4.2 Характеристики измерения

ΠΑΡΑΜΕΤΡ	ЗНАЧЕНИЕ						
Время измерения: одноволновое измерение двухволновое измерение	< 15 c < 20 c						
Диапазон длин волн: стандартный	400 - 750 нм						
Диапазон измерения: 400 - 750 нм	0 - 4,000 OD						
Разрешение экрана:	0,0001 OD						
Точность: 450, 492 нм 0,000 - 2,000 OD 2,000 - 3,000 OD	≤ (1.0 % + 0.010 OD)* ≤ (1.5 % + 0.010 OD)*						
Сходимость: 450, 492 нм 0,000 - 2,000 OD 2,000 - 3,000 OD	≤ (0.5 % + 0.005 OD)* ≤ (1.0 % + 0.005 OD)*						
Линейность: 450, 492 нм 0,000 - 2,000 OD 2,000 - 3,000 OD	≤ 1 % ≤ 1.5 %						
Выбор длины волны: стандартный фильтр	Узкополосные интерференционные фильтры. На диск фильтров можно установить до 8 фильтров.						
Точность длины волны фильтра:	Центральная длина волны ± 2 нм						
Ширина полосы фильтра: при 50%-ном пропускании	10 ± 2 нм						
Источник света:	Светодиод						
Компьютерный интерфейс:	USB						
Все подключенные устройства должны соответствовать и входить в перечень согласно нормативам IEC 60950-1 «Оборудование для							

информационных технологий — безопасность» и аналогичным местным стандартам.

\* лучше или равно сумме x % от измеренного значения и значения OD с соответствующим коэффициентом



### 4.4.3 Микропланшеты

В фотометре INFINITE F50 PLUS могут применяться все 96-луночные микропланшеты с прозрачным дном (плоские, С-, U- и V-образные; включая лунки в стрипах), если они отвечают следующим стандартам:

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



#### ПРЕДОСТЕРЕЖЕНИЕ ИСПОЛЬЗУЙТЕ ТОЛЬКО МИКРОПЛАНШЕТЫ БЕЗ КРЫШЕК. НЕ ИСПОЛЬЗУЙТЕ МИКРОПЛАНШЕТЫ ВЫСОТОЙ БОЛЕЕ 15,2 ММ.

#### Работа с микропланшетом

Вставлять или извлекать микропланшет разрешается только при полностью извлеченном транспорте планшета (как показано ниже) и остановленном электродвигателе перемещения транспорта. Не открывайте крышку корпуса, если индикатор состояния горит красным светом.



#### ПРЕДУПРЕЖДЕНИЕ ДЛЯ РАБОТЫ С МИКРОПЛАНШЕТОМ ВСЕГДА НАДЕВАЙТЕ ЗАЩИТНУЮ ОДЕЖДУ И ОДНОРАЗОВЫЕ ПЕРЧАТКИ.



Положение лунки А1 Транспорт планшета – полностью извлечен

## 4.5 Принадлежности прибора

В приведенном ниже списке представлены дополнительные принадлежности фотометра INFINITE F50 PLUS:

- дополнительные фильтры;
- инструмент для монтажа фильтров;
- планшет MultiCheck<sup>™</sup> для приборов семейства INFINITE F50

Для получения дополнительной информации и проверки возможности поставки интересующих принадлежностей обратитесь к местному торговому представителю Tecan.



## 5. Контроль качества

## 5.1 Введение



#### ПРЕДОСТЕРЕЖЕНИЕ В СЛУЧАЕ ВОЗНИКНОВЕНИЯ СОМНЕНИЙ В АНАЛИТИЧЕСКИХ ХАРАКТЕРИСТИКАХ INFINITE F50 PLUS ПОЛЬЗОВАТЕЛЬ ДОЛЖЕН СЛЕДОВАТЬ ИНСТРУКЦИЯМ ПО КОНТРОЛЮ КАЧЕСТВА ИЛИ ОБРАТИТЬСЯ В МЕСТНЫЙ СЕРВИСНЫЙ ЦЕНТР ТЕСАN.

В этой главе приводятся сведения о процедуре самопроверки прибора и инструкции по простой проверке качества его работы.

## 5.2 Процедура самопроверки

При подключении фотометра INFINITE F50 PLUS к управляющему программному обеспечению Magellan выполняется проверка электродвигателей и датчиков с инициализацией транспорта планшета и диска фильтров.

Перед каждым измерением проводится калибровочная процедура самопроверки, предназначенная для проверки надлежащей работы прибора и выполнения калибровки оптической системы.

## 5.3 Эксплуатационная квалификация (ЭК)

Проверить надлежащую работу прибора и точность получаемых результатов можно с помощью приведенных ниже тестов.

Воспроизводимость и точность показаний прибора может меняться в зависимости от исследуемого раствора и используемого микропланшета.

Для предотвращения этого эффекта на заводе все приборы тестируются при помощи калибровочного планшета, что позволяет устранить влияние характера исследуемого раствора и размещения микропланшета на точность результатов исследований.

## 5.3.1 Tecm MultiCheck

Тест MultiCheck представляет собой автоматизированную проверку характеристик прибора, включая проверку точности, линейности, сходимости измерений и соответствия стандартам, контролепригодным согласно требованиям NIST (Национального института стандартов и технологий США).

### 5.3.2 Тестирование микропланшета

Если лунки микропланшета имеют различную оптическую плотность, это повлияет на результаты измерений.

Проверить неравномерность оптической плотности ячеек можно посредством измерения на пустом микропланшете.

Значения оптической плотности, полученные на пустом микропланшете, должны колебаться в узком диапазоне. Например: ± 0,010 OD.



Если значения оптической плотности выходят за указанные пределы, использовать микропланшет нельзя.

При использовании двухволновых измерений влияние отклонений оптической плотности лунок на результаты устраняется либо уменьшается до приемлемых пределов.

#### 5.3.3 Тест сходимости прибора на образцах жидкости

Эта процедура используется для проверки сходимости измерений. Рекомендуется использовать микропланшет с плоским дном.

Заполните новый микропланшет только что приготовленным раствором оранжевого Ж с разным разведением для разных лунок, чтобы значения оптической плотности получились в широком диапазоне. В лунках должно быть не менее 200 мкл раствора. Серии разведения должны соответствовать значениям оптической плотности от 0,1 до 3,0 OD. Значение оптической плотности 3 OD получается при концентрации оранжевого Ж (номер по каталогу компании Sigma: O7252) приблизительно 125 мг·л<sup>-1</sup>.

Задайте тестовый прогон с использованием фильтра 492 нм и измерьте микропланшет не менее трех раз.

Для каждой лунки вычислите следующие параметры:

- среднее значение оптической плотности;
- стандартное отклонение.

#### Примеры

#### Показания от 0,000 до 2,000 OD

Стандартное отклонение для каждой лунки должно находиться в пределах (0,5 % + 0,005 OD).

Расчет максимально допустимого отклонения для среднего значения оптической плотности 1,000 OD:

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

#### Показания от 2,001 до 3,000 OD

Стандартное отклонение для каждой лунки должно находиться в пределах (1,0 % + 0,005 OD).

Расчет максимально допустимого отклонения для среднего значения оптической плотности 2,400 OD:

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

#### Показания выше 3,000 OD

Показания выше 3,000 OD используются только в ориентировочных целях. Гарантировать сходимость таких значений невозможно.



## 5.3.4 Тест линейности прибора на образцах жидкости

Линейность измерений прибора и программного обеспечения на используемой длине волны можно проверить с помощью серии разведений раствора.

Результат теста зависит от чистоты использованного красителя и величины мениска жидкости в лунках.

В качестве контрольного раствора для измерений на длине волны 492 нм можно использовать серию разведений раствора оранжевого Ж.

Серии разведения должны соответствовать значениям оптической плотности от 0,1 до 3,0 OD. Значение оптической плотности 3 получается при концентрации оранжевого Ж (номер по каталогу компании Sigma: O7252) приблизительно 125 мг·л<sup>-1</sup>.

Для проверки на других длинах волн следует использовать другие растворы.

В лунки микропланшета вносят 200 мкл каждого раствора, для устранения погрешности, связанной с неточностью дозировки, следует внести раствор каждой концентрации не менее чем в две лунки.

Затем выполняется измерение и строится линейная регрессия средних значений оптической плотности в зависимости от концентрации раствора.

Определите невзвешенный коэффициент детерминации R<sup>2</sup> линии регрессии.

Обычно невзвешенный коэффициент детерминации для стандартного применения равен или лучше R<sup>2</sup> = 0,998.



Примечание Возможны отклонения полученных значений вследствие погрешности дозировки раствора.



## 6.1 Введение

Файлы примера, основанные на методах и рабочих пространствах Magellan, предназначены для ознакомления с этим программным обеспечением и облегчения работы с ним. Файлы примера количественного и качественного анализа ELISA устанавливаются автоматически при установке Magellan.

# 6.2 Пример с пошаговыми инструкциями: количественный анализ ELISA

В этой главе рассмотрен пример создания метода количественного анализа с пошаговыми инструкциями. Следуя инструкциям из описания диагностического комплекта, вы научитесь определять оценки.



Примечание В Magellan файлы примера автоматически отображаются в списке методов. В Magellan Tracker эти файлы находятся в задаваемом по умолчанию каталоге и должны быть конвертированы.

## 6.2.1 Описание диагностического комплекта

В предоставляемом производителем описании диагностического комплекта, в котором рассматриваются вопросы количественного анализа иммуноглобулина М, обнаружения антител и анализа ELISA, приведены следующие инструкции: Схема планшета

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

BLK = бланк, NC = отрицательный контроль, C1 - C6 = стандарты (калибраторы),

S1 - S... = образцы



#### Измерение и оценка

Измерьте планшет, используя длину волны 492 нм и опорную длину волны 620 нм.

В позиции лунки А1 используется бланк.

Концентрации стандартов (калибраторов):

Стандарт 1	5 УЕ/мл
Стандарт 2	10 УЕ/мл
Стандарт 3	20 УЕ/мл
Стандарт 4	40 УЕ/мл
Стандарт 5	80 УЕ/мл
Стандарт 6	160 УЕ/мл

После учета начального уровня строится график зависимости оптической плотности (492 - 620 OD) от концентрации. Проходящая через полученные точки линия регрессии представляет собой калибровочную кривую.

Интерпретация результатов измерения:

IgM < 18 УЕ/мл	Отрицательный
18 УЕ/мл ≤ IgM < 22 УЕ/мл	Промежуточный
IgM ≥ 22 УЕ/мл	Положительный

Вычисленная концентрация иммуноглобулина М (IgM) обоих отрицательных контролей должна быть менее 8 УЕ/мл.

#### Обработка данных

По завершении измерения автоматически сохраняется файл данных (рабочее пространство) и создается отчет, в котором отражаются параметры измерения, схема планшета, значения для учета нулевого уровня, калибровочная кривая, концентрации иммуноглобулина М, граничные значения, качественные результаты измерения образцов и результаты проверки.

Кроме того, схема планшета и качественные результаты сохраняются также в виде файла ASCII.



## 6.2.2 Создание метода

В диалоговом окне Список мастеров выберите Созд./редактир. метода и нажмите ОК. На странице Добро пожаловать! мастера Создание/редактирование метода нажмите кнопку Продолжить для открытия диалогового окна Выбор файла. Выберите Создать.

Создание/редактирование метода				×
<ul> <li>Создать</li> <li>Открыть</li> </ul>	Показать:	Файлы с данного прибора 💌	Предварит. просмотр	
🗁 mth	Name 🔺	Комментарии	Состояние	
Справка Отмена <<< Назад		Сделайте Е	Заш выбор	
		©	2009 Tecan	

#### Параметры измерения

Нажмите кнопку Сделайте Ваш выбор для открытия страницы Параметры измерения.





На панели **Длина волны** выберите в поле Длина волны измерения значение 492 нм, а в поле Референсная длина волны — значение 620 нм.

🍦 🔻 Оптич	еская плотнос	ть		3
_Длина волны			Метка	
Измерение:	492 nm	~	Имя: Метка1 💌	
🗹 Эталон:	620 nm	*		

Для продолжения работы мастера нажмите кнопку **Выбрать параметры** измерения. Отображается окно **Схема планшета**.

Создание/редактирование метода														×
Отмена Поетор Выбрать все неисполь	3 10% 1	100% <b>1</b> 98	елич											
Схена нетода 🛞														
ть/редакти	0	1	2	3	4	5	6	7	8	9	10	11	12	
ровать мето	A											Назначение л	Эксп. группа	
A	B											ST PC	Идент. №.	
	C											NC LPC HPC	Повторения	ור
	D											RF 🗸	Bce	D:
	E											Выбор загрузки	Удаление загрузки	
	F													
	G													AN Austria
	H													TEC
Справка Отмена <<< Назад													Далее	



#### Определение схемы

Задайте схему планшета в диалоговом окне Назначение лунки, расположенном на правой стороне экрана.

Назн	Назначение лунки								
сид	дентиф	ик. —	Эксп. группа:						
	SM	^	1						
	BL		_Идент. №.						
	ST		≠ 4						
	PC								
	NC								
	LPC		Повторения						
	HPC		≠ 4						
	BF		💿 Зак. номер:						
	RF	~	1						
	ред. и,	дент.	OBce						
	Выбор загруз	р ки	Удаление загрузки						

В группе Идентификатор выберите значение BL (бланк).

В группе Эксп. группа остается число 1.

В группе Повторения автоматически включается радиокнопка Все.

Щелкните на лунке А1. Рамка лунки станет красной.

Нажмите кнопку Заполнить выбранную область . Лунка помечается идентификатором выбранного типа.



#### Примечание

Отдельную лунку можно также заполнить двойным щелчком на ней.

Теперь выберите в диалоговом окне **Назначение лунки** следующие параметры:

В группе **Идентификатор** выберите значение **NC** (отрицательный контроль).

В группе Эксп. группа остается число 1.

В группе Повторения автоматически включается радиокнопка Все.

Щелкните на лунке **B1** и, не отпуская кнопки мыши, переведите указатель к лунке **C1**. Теперь красной рамкой обведены лунки **B1** — **C1**.

Нажмите кнопку Заполнить выбранную область. Лунки помечаются идентификатором выбранного типа.

Определим лунки **D1** — **G2** под заполнение стандартами (калибраторами). Теперь выберите в диалоговом окне **Назначение лунки** следующие параметры:

В группе Идентификатор выберите ST (стандарт).



В группе Эксп. группа остается число 1.

В группе Повторения выберите между Фикс. число и Все.

Фикс. число:

Этот параметр доступен только для стандартов и образцов, на которых можно использовать идентификаторы.

Если радиокнопка **Фикс. число** включена, в ее поле можно ввести число. Данное число определяет, сколько повторений будет иметь данный метод. Введенное число повторений действует в выбранных лунках в отношении всех идентификаторов. Следовательно, число выбранных лунок должно быть кратно введенному числу повторений.

#### Bce:

Все выбранные лунки задаются как повторения. Если для выделенных лунок выбран существующий идент. номер, они будут добавлены в качестве повторений к уже существующим повторениям. Выделенные лунки с любыми другими типами идентификаторов также будут добавлены в качестве повторений к уже существующим повторениям.

Две кнопки со стрелками задают направление нумерации повторений (горизонтальное или вертикальное).

Включим для этого примера радиокнопку Фикс. число и зададим число 2.

В окне Идент. №. и группе Повторения выберите вертикальные стрелки.

Затем выберите лунки D1 — G2 и нажмите кнопку Заполнить выбранную область.



Примечание Выбор лунок выполняется следующим образом: щелкните на лунке D1 и, не отпуская кнопки мыши, переведите указатель к H1. Затем, удерживая нажатой клавишу Ctrl, перетащите

указатель от А2 к G2.



Создание/редактирование мет	года														X
Отмена Повтор Выбрать все неи	спольз.	10%	100% 🗌 9a	елич											
Скена нетода 📀	(e)														
Конц, Разв, Эталонз	арть/редакти	0	1	2	3	4	5	6	7	8	9	10	11	12	$\sum$
Концентрации 🛞	ровать мето	A	BL1 1/1	ST1_3 2/2									Назначение л Идентифик. SM	унки Эксп. груг 1	ina:
Оценивать данные 🔗 Определение граничных Проверка контроля качес	A	B	NC1 1/2	ST1_4 1/2									ST PC	Идент. №	
<ul> <li>Обработка данных (※)</li> <li>Экспорт данных</li> <li>Печатный отчет</li> </ul>		C	NC1 2/2	ST1_4 2/2									LPC HPC	Повторен	
Автоматическая обработ     Дополнительно     Запросы пользователя		D	ST1_1 1/2	ST1_5 1/2									RF 🗸	© Зак. ног 2 	10p)
Формат чисел Принечания к нетоду		E	ST1_1 2/2	ST1_5 2/2									Выбор загрузки	Удалені загрузк	ле И
		F	ST1_2 1/2	ST1_6 1/2	,										
		G	ST1_2 2/2	ST1_6 2/2											AN Austria
		H	ST1_3 1/2												TEO
	Į														
		_	_	_		_	_	_	_	_	_			_	_
Справка Отмена <<< Назад														Далее	

Теперь окно Схема планшета выглядит следующим образом:

Щелкните на панели инструментов кнопку **Выбрать все** неиспользованные для выбора всех пустых лунок шаблона. Затем, удерживая нажатой клавишу Ctrl, щелкните на лунке **H12**, чтобы она осталась пустой и непомеченной.

В группе Идентификатор диалогового окна Назначение лунки выберите SM (образец).

В группе Эксп. группа остается число 1.

В группе Повторения выберите Фикс. число и 2.

В окне **Идент. №.** и оставьте значение 1, а в группе **Повторения** выберите **вертикальные стрелки**. Нажмите кнопку **Заполнить выбранную область**. Процедура определения схемы планшета завершена.

#### Преобразования

Выберите в группе **Преобразованные данные** панели управления на левой стороне окна пункт **Добавить новое преобразование...**, чтобы определить учет нулевого уровня.



Отображается диалоговое окно с запросом о необходимости определения учета нулевого уровня. Нажмите кнопку **Да**. Отображается следующее окно:

Создание/редактирование метода	í.														
Отмена Повтор	10		100% 🗌 Ув	елич											
Ś	Вход.		Разностные да	нные			~						1	Константы	Паранетры
Пурма мето ва		fx	x-BL1							~	Инеющиеся	данные (неско	лько наб 🔻	Функции и К	онстанты 🔻
🕉 Схема планшета															
Конц, Разв, Эталонз															
Чо Преобразованные 🔊 🔮															
Учет нулевого уровня	6		1	2	3	A	5	ß	7	8	0	10	11	12	
Добавить новую трансфо							0		u			10			
🕎 Концентрации 🛞 🖁			BL1	ST1 3	SM1 1	SM1 5	SM1 9	SM1 13	SM1 17	SM1 21	SM1 25	SM1 29	SM1 33	SM1 37	
Калибровочная кривая			1/1		2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	
Оценирать данные			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Определение граничных			NC1	ST1_4	SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
Проверка контроля качес			1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	1/2	
/ Обработка данных 🛞			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
Экспорт данных	0	-	NC1	ST1_4	SM1_2	SM1_6	SM1_10	SM1_14	SM1_18	SM1_22	SM1_26	SM1_30	SM1_34	SM1_38	
Печатный отчет	0	9	2/2 x BL1	2/2 v RI 1	2/2 x BL1	2/2 x BL1	2/2 x BI 1	2/2 × BL1	2/2 x RI 1	2/2 x BL1	2/2 x BL1	2/2 x BL1	2/2 x BL1	2/2 x BL1	
Автопатическая оорасот			CT4 4	OTA E	Chill O	Child 7	Child dd	Child dE	Child do	Child 00	Ch44 07	Child Od	Child DE	Child 20	
💥 Дополнительно 🛞	l f		1/2	511_5 1/2	SMIL_3	SMT_/	5MI_II 1/2	5MI1_15 1/2	5MI_19 1/2	5NII_23	5WI1_27	5WI_3T	5WI1_35 1/2	5MT_39 1/2	
Запросы пользователя		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	
Примечания к методу			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	
		_	ST1_2	ST1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	
		F	1/2		1/2	1/2	1/2		1/2	1/2	1/2		1/2	1/2	
			x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	x-BL1	
			ST1_2	ST1_6	SM1_4	SM1_8	SM1_12	SM1_16	SM1_20	SM1_24	SM1_28	SM1_32	SM1_36	SM1_40	Strig
	0	5)	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	2/2	1 At
			X-BLI	X-BLI	X-BLI	X-BLI	X-BLI	X-BLI	X-BLT	X-BLI	X-BLT	X-BLI	X-BLI	X-DL1	CAP
	I B	1	SI1_3 1/2	SM1_1	SM1_5	SM1_9	SM1_13	SM1_17	SM1_21	SM1_25	SM1_29	SM1_33	SM1_37		Ψ
	0	U	x-BI 1	x-BI 1	x-BL1	x-BL1	x-BI 1	x-BI 1	x-BI 1	x-BI 1	x-BI 1	x-BI 1	x-BI 1		
			_	_	_	_	_	_		_	_	_	_		
Справка															
Отмена <<< Назад														Далее	
													(D. 200	19 Teran	

В поле Входные данные автоматически выбирается значение Разностные данные Если ранее был подтвержден выбор определения учета нулевого уровня, программа автоматически задает имя Учет нулевого уровня в качестве названия преобразования на панели управления.

В окне **Формула** автоматически появляется значение **x-BL1** для данного учета нулевого уровня, где «х» означает текущее входное значение лунки, а «BL1» — среднее по экспериментальной группе 1 значение для лунок с бланками.

Для получения дополнительной информации об определении и назначении преобразований см. Руководство пользователя Magellan.

В каждой лунке отображается некоторая информация, значение которой, рассматриваемое на примере лунки А5, приведено в таблице ниже.

SM1_9	Образец, экспериментальная группа 1, номер идентификатора образца 9.
2/2	Номер повторения – 2, общее число повторений – 2.
x-BL1 или 1	Назначено преобразование x-BL1 (если выделено «Преобразование») или Степень разведения 1 (если выделено «Конц, Разв, Эталонзначения»).





#### Определение концентрации, разведения и эталонного значения

Выберите в группе **Схема метода** на панели управления пункт **Конц.-, Разв.-, Эталон.-значения**, чтобы задать значения согласно описанию диагностического комплекта.

Стандарт 1	5 УЕ/мл
Стандарт 2	10 УЕ/мл
Стандарт 3	20 УЕ/мл
Стандарт 4	40 УЕ/мл
Стандарт 5	80 УЕ/мл
Стандарт 6	160 УЕ/мл

Убедитесь, что в раскрывающемся списке **Выбор идентификатора** отображается значение **ST**.

В столбце **Идентификатор** отображается список стандартов из экспериментальной группы 1. Введите в столбце **Концентрация** строки **ST1\_1** число **5**, а в поле **Единица** — «УЕ/мл». Введите в столбце **Концентрация** строки **ST1\_2** число **10**. Единица измерения определяется только один раз и распространяется на все стандарты. Аналогичным образом введите значения для стандартов ST1\_3 — ST1\_6.

Отображается экран со схемой планшета и концентраций:





#### Калибровочная кривая

Для задания калибровочной кривой выберите в группе Концентрации пункт Калибровочная кривая.

Приведенная ниже информация содержится в описании диагностического комплекта.

После учета начального уровня строится график зависимости оптической плотности (492 - 620 OD) от концентрации. Проходящая через полученные точки линия регрессии представляет собой калибровочную кривую.

На вкладке **Данные** выберите в расширяющемся списке «Входные данные» значение **Учет нулевого уровня**.

анные Тип анализа Отрезки Ось График
Вход. данные: Учет нулевого уровня
<ul> <li>Стандарты из схемы</li> </ul>
О Стандарты из внеш. файла: Выбрать
🔿 Стандраты из экпериментальной группы: 🔍
🔾 Без графика калибровочной кривой
Дополнительные концентрации

На вкладке Тип анализа выберите значение «Линейная регрессия».

<ul> <li>Линейная регрессия</li> <li>Нелинейная регрессия</li> </ul>	Масштаб, данных: Lin(x)Lin(y) 🗸
○ Кубический сплайн ○ <u>А</u> kima	and Amount 2
О Полиномиальный 2 🗘	
○ Четыре параметра	
O Marquardt с четырьмя параметрами	Acrony
○ 5 Параметров	donom
○ LogitLog	
1	

2023-06



Метка: Концентрация [UA/n	n[]
Цвет:	🗌 Логарифм. масштаб
💿 Автомат. выбор диапазона	
🔿 Диапазон	Мин.: Макс.:
🗹 Сетка	Цвет:
Ось Ү	
Метка: Учет нулевого уров	ня
Цвет:	🗌 Логарифм. масштаб
<ul> <li>Автомат. выбор диапазона</li> </ul>	
О диапазон	Мина: Макса
_	

На вкладке Ось определите маркировку и масштабирование оси.

На вкладке **График** определите название графика, параметры кривой, шрифт и отображение различных параметров графика.

Метка: IgM-ELISA			
Цвет:	T		
Кривые			
	Метка: Гр. 1		
Цвет:	Скрыть кривую		
Символ:	Толщина линии: 🛛 😂		
Шрифт	Показать		
💿 Малый	🗹 Легенда	🗹 Отрезки	
<ul> <li>Средний</li> <li>Большой</li> </ul>	🗹 Базовые точки	🗹 Планки погрешностей	



#### Определение граничных значений

Для задания границ качественной оценки выберите в группе **Оценка данных** на панели управления пункт **Определение граничных значений**.

В описании диагностического комплекта приведены следующие инструкции:

Интерпретация результатов измерения:

IgM < 18 УЕ/мл	Отрицательный
18 УЕ/мл ≤ IgM < 22 УЕ/мл	Промежуточный
IgM ≥ 22 УЕ/мл	Положительный

Определение граничных условий производится следующим образом:

В списке «Входные данные» выберите значение Средняя конц. (УЕ/мл).

Таблица **Граничные значения** представляет собой шкалу, на которой указаны верхнее и нижнее пороговые значения для столбцов **Пределы** и **Метки**. В столбце **Пределы** введите 22 в качестве первого (верхнего) предела и 18 — в качестве второго (нижнего) предела.

Введите в отдельные ячейки столбца **Метки** интерпретации результатов (**Положительный**, **Промежуточный** и **Отрицательный**). Выберите из раскрывающейся палитры следующие цвета:

- Положительный красный
- Промежуточный синий
- Отрицательный зеленый

Отображается следующий экран:

Вход. данные: Средняя конц. (UA/ml)	
Граничные значения	
Цвета Метки Пределы positive 22 intermediz negative 18 regative 4 intermediz interm	
Веод формулы Переменная Операторы Функции BL1 • + • and •	Конкурентный тест Конкурентный тест Зыбор результ. примен. грани. значений

Для выбора типов идентификатора, для которых требуется показать результаты применения граничных условий, нажмите кнопку **Выбор результ. примен. грани. значений**.



#### Определение проверок контроля качества

Выберите в группе **Оценка данных** на панели управления пункт **Проверки** контроля качества. Для гарантии достоверности результатов тестирования необходимо определить критерии проверки.

В этом примере должно быть выполнено следующее требование:

Вычисленная концентрация иммуноглобулина М (IgM) обоих отрицательных контролей должна быть менее 8 УЕ/мл.

В списке Входные данные выберите значение Единичн. конц. (УЕ/мл).

В первой строке введите NC1\_1<8



Примечание «NC1\_1» означает «отрицательный контроль экспериментальной группы 1, повторение 1».

Во второй строке введите NC1\_2<8.

Теперь диалоговое окно **Проверки контроля качества** отображается в следующем виде:

сдини				
Условия проверки				
NC1_1<8				
NC1_2<8				
од формулы				
Попомоннов				
переменная	Операторы	Функции		
BL1	Операторы	о Функции	~	
BL1	Операторы	Функции and	~	
ВL1	Операторы + паншета к планшету	о Функции	<b>v</b>	
ВL1	Операторы +	о Функции		
ВL1 всодные данные:	Операторы + паншета к планшету Разностные данны	Функции and	<ul> <li></li> <li></li> </ul>	
ВL1	Операторы + + паншета к планшету Разностные данны	е Функции	<ul> <li></li> </ul>	 
ВL1 всодные данные: Управление 1:	Операторы + паншета к планшету Разностные данны ВL1 Сре	е еднее значение:	<b>×</b>	Ť
ВL1 ВL1 ВСЛИВНИ ВСЛИВИИ ВСЛИВИ ВСЛИВИ ВСЛИВИ ВСЛИВИ ВС ВСЛИВИВИ ВСЛИВИ ВСЛИВИ ВСЛИВИ ВСЛИВИ ВСЕЛИВИ ВСЕЛИВИИ ВСЛИВИИ ВСЛИВИИ ВСЛИВИИ ВСЕЛИВИВИ ВСЕЛИВИВИ ВСИВИИ ВСЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИВИ ВСИВИИ ВСЕЛИВИИ ВСИВИ ВСИВИИ ВСЕЛИВИ ВСЕЛИВИ ВСЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЛИВИ ВСЕЛИВИ ВСЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЛИВИИ ВСЛИВИ ВСЛИВИ ВСЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЕЛИВИ ВСЛИВИ ВСТИВИ ВСЕЛИВИ ВСЕЛИВИ ВСТИ ВСИВИ ВСЛИВИ ВСЛИВИ ВСТИВИ ВСТИВИ	Операторы + паншета к планшету Разностные данны ВL1 Сре ВL1 Сре	Функции Функции аnd е е е е е е е е е е е е е е е е е е е	••••••••••••••••••••••••••••••••••••••	Ť
ВL1 ВL1 ВСЛИВНИКАЧЕСТВА ОТ П Входные данные: Управление 1: Управление 2:	Операторы + наншета к планшету Разностные данны ВL1 Сре ВL1 Сре ВL1 Сре	Функции Функции and e e e e e e e e e e e e e e e e e e e		Ť
ВL1 ВL1 ВССТВА ОТ П Входные данные: Управление 1: Управление 2: Управление 3:	Операторы + наншета к планшету Разностные данны ВL1 Сре ВL1 Сре ВL1 Сре	Функции аnd е е еднее значение: еднее значение: еднее значение:	CO:	Î.
ВL1 ВL1 Всодные данные: Управление 1: Управление 2: Управление 3: Управление 4:	Операторы 	Функции аnd e е е е е е е е е е е е е е	CO: CO: CO: CO:	1
ВL1 ВL1 ВL1 ВСС СТВА ОТ П Входные данные: Управление 1: Управление 2: Управление 3: Управление 4: О Все рабочие про	Операторы 	Функции аnd е е е е е е е е е е е е е	CO: CO: CO: CO:	<u>↑</u>
ВL1 ВL1 ВL1 ВСОДНЫЕ ДАННЫЕ: Управление 1: Управление 2: Управление 3: Управление 4: О Все рабочие про Рабочее простое	Операторы аншета к планшету Разностные данны ВL1 Сре ВL1 Сре ВL1 Сре ВL1 Сре ВL1 Сре ВL1 Сре ВL1 Сре ВL1 Сре	• Функции • алd • алд •	 CO: CO: CO: CO:	<u>↑</u>



### 6.2.3 Настройка печатного отчета

Выберите в группе **Обработка данных** на панели управления пункт **Печатный отчет**. Отображается следующий экран:

Печатать н	как писок ОМатрица		Вы можете Выберите од список выбра	использовать функцию drag_drop, чтобы со ин пункт из имеющихся данных в списке и п нных данных. Поместите его в соответству отчете.	оздать отчет. ередвинте его в ющую позицию в
Доступные д	анные:			Выбранные данные:	
	нные прибора еньшенные данные Разностные данные Разностные данные - Разностные данные - еобразованные данные - еобразованные данные нцентрации личественные результа сент. образцов ема метода итерии проверки контр раметры измерения мментарии ирал ошибок ирал ошибок длись	Сред Стан, Коэф аты оля к.	Вставить Добавить	<ul> <li>Е. Сбор данных</li> <li>Параметры измерения</li> <li>Вертик, список</li> <li>Вертик, список</li> <li>Разностные данные</li> </ul>	Вверх Вниз Свойства Удалить Удалить Все

Все доступные для отчета данные содержатся в окне **Доступные данные** на вкладке **Выбор данных**. Для переноса данных в окно **Выбранные данные** используются кнопки **Вставить и Добавить**. Это можно также сделать перетаскиванием.

На панели **Печатать как** выберите между печатью данных в виде матрицы или в виде списка с заданной ориентацией.

В этом примере создадим отчет, содержащий параметры измерения, схему планшета, значения для учета нулевого уровня, калибровочную кривую, концентрации иммуноглобулина М, граничные значения, качественные результаты измерения образцов и результаты проверки.

Перед созданием отчета из окна Выбранные данные необходимо удалить заданный по умолчанию параметр Вертик. список/Разностные данные. В окне Выбранные данные остается только элемент Параметры измерения. На панели «Печатать как» вместо радиокнопки Список необходимо включить радиокнопку Матрица.



Печатать как О Список 💿 Матрица	ы ножете использовать функцино drag drop, чтобы создать отчет. Выберите один пункт из имеющихся данных в списке и передвинте его в список выбранных данных. Поместите его в соответствующую позицию в отчете.
Доступные данные:	Выбранные данные:
Даняве просра     Хуменьшенные данные     Разностные данные - Сред     Разностные данные - Стан,     Разностные данные - Стан,     Разностные данные - Стан,     Разностные данные - Кооф     Концентрации     Количественные результаты     ', Идент, образцое     Схема иетода     Критерии проверки контроля к.	Ставить В 2 Сор даннол Добавить Вно Свойства. Удалить во
Параметры измерения Комментарии Журнал ошибок Журнал контроля Подпись	Инпорт

В окне **Доступные данные** выберите элемент **Схема метода/Схема** и прикрепите его к отчету в виде матрицы нажатием кнопки **Добавить**. Затем вставьте в матрицу элементы **Учет нулевого уровня**, **Средняя конц.** (УЕ/мл) и **Результаты применения граничных значений**, для чего необходимо выбрать их и нажать кнопку **Вставить**.

Добавьте к выбранным данным элементы График: Калибровочная кривая, Определение граничных значений и Критерии контроля качества. Теперь этап настройки данных в процедуре определения отчета завершен; диалоговое окно Печатный отчет выглядит следующим образом:



На вкладках **Верхний колонтитул** и **Нижний колонтитул** определите схему верхнего и нижнего колонтитулов отчета (подробнее см. Руководство пользователя Magellan).

#### Экспорт данных

Выберите в группе **Обработка данных** на панели управления пункт **Экспорт данных**. В этом примере сохраним схему и результаты применения граничных значений в виде файла ASCII. Выберите элементы **Схема** и **Результаты применения граничных значений** в окне **Доступные** 



**данные** и нажмите стрелку →, чтобы вставить их в окно **Выбранные данные**. Отображается следующий экран:

🕀 🛄 Данные прибора	^		Схема	Вверх
🖃 🤆 Уменьшенные данные			Разностные данные	Durin
				БНИЗ
Разностные данные - Среднее зн				
📃 Разностные данные - Стандартн				
📳 Разностные данные - Коэффициє				
🗄 🔏 Преобразованные данные				
🗉 🚺 Концентрации		->		
Результаты применения граничных за				
н <sup>1</sup> 2, Идент. образцов		<-		
🖃 📶 Схема метода				
Позиции лунок				
📃 Имена метода для стрипов	-			
📳 Первоначальные концентрации				
	-			
				1



Примечание Элементы «Схема» или «Список идентификаторов образцов» должны всегда включаться в экспортируемые данные.

#### Автоматическая обработка данных

Выберите в группе **Обработка данных** на панели управления пункт **Автоматическая обработка данных**.

Автоматически	
🔲 Загрузить список ID образцов	Дополн
✓ Сохранить рабочее пространство	Дополн
Экспорт в ASCII файл	
🗌 Экспорт в ASTM файл (LIS)	Дополн
Экспорт в Excel	
Печать	
🗹 Просматривать результ. после измерения	Дополн

Выберите элементы Экспорт в ASCII файл и Просматривать результ. после измерения. В Magellan Tracker по умолчанию выделен элемент Сохранить рабочее пространство, который не может быть изменен.



#### Сохранение метода

Нажмите **Далее**. Откроется окно **Сохранить как**. Введите имя файла метода и при необходимости заполните остальные поля.

Создание/редактирование метода			
Папка: 🗀 mth	Name 🔺	Комментарии	Состояние
Имя файла: Метод1.mth			
Замечание к файлу:			
Комментарии для контрольного журнала:			
		_	
Упоряд. избранное Журнал контроля	Подписи	Пароль метода:	
Справка Отмена (<< Назад Зап	устить этот метод сейчас	Сохранить и законч	ять 🕨
		© 2009 T	ecan

<b>Имя файла</b> текстовое поле	Обязательный параметр. Предлагаемое системой имя файла можно изменить.
Комментарии к файлу текстовое поле	Введенные здесь комментарии сохраняются и отображаются вместе с именем файла.
Комментарии для контрольного журнала текстовое поле	Введенные здесь комментарии сохраняются в контрольном журнале. Данный параметр доступен только в <b>Magellan Tracker</b> .
Упорядочить избранное кнопка	Открытие диалогового окна <b>Упорядочить</b> избранное.
Пароль метода текстовое поле	Введите пароль для защиты метода.
Запустить этот метод сейчас флажок	Немедленный запуск метода после нажатия кнопки «Сохранить и закончить».

Подробнее см. Руководство пользователя Magellan.



## 6.2.4 Запуск метода

Если в диалоговом окне **Сохранить как** мастера **Создание/редактирование метода** установлен флажок **Запустить этот метод сейчас**, после нажатия кнопки **Сохранить** отображается мастер/диалоговое окно **Запуск измерения**. В диалоговом окне **Запуск измерения** отображается задаваемое по умолчанию имя рабочего пространства, которое можно изменить.

Запуск измерения		Σ
Измерение Рабоч. простр: 06112009-001.wsp Метод: Метод1.mth Список идент. образцое Испол-зовать податчик Загрузить планшет Регул. температ Перемещения Контр. впрыска	Кинет. с произв. инт Изменить схему Вставить Текущие: Не ог °С Цель: Не ог °С Настр. полож. по верт	Параметры измерения Планшет Описание планшета: [GRE96ft] - Greiner 96 Fl. Планшет с крышкой: Нет Штриккоа: Нет Часть планшета Диапазон: A1:H12 Оттическая плотность Длина волны измерения: 492 nm Опорная длина волны: 620 nm Метка: Метка1
Справка		Пуск
		@ 2009 Tecan

Для запуска измерения нажмите **Пуск**. Рабочее пространство создается автоматически и содержит все ранее введенные сведения. В нем сохраняются все значения измерений. Ход измерения отражается в диалоговом окне состояния.

По завершении измерения открывается диалоговое окно **Результаты**, в котором можно просмотреть все результаты и расчеты. При выполнении измерения без необходимых жидкостей (например, стандартов) могут появиться сообщения об ошибках.



### 6.2.5 Оценка результатов

Для просмотра и оценки исходных данных выберите элемент **Оценка результатов**. Также возможны просмотр параметров оценки данных и повторная оценка данных.

В этом разделе приведено описание мастера **Оценка результатов** на примере демонстрационного файла рабочего пространства, устанавливаемого одновременно с Magellan.

В диалоговом окне *Список мастеров* выберите мастер **Оценка результатов**.

На странице **Добро пожаловать!** мастера **Оценка результатов** нажмите кнопку **Далее** для открытия диалогового окна **Выбор файла**.

Выберите из списка файлов рабочее пространство Quantitative ELISA example\_Sunrise\_InfiniteF50.wsp и нажмите кнопку Сделайте Ваш выбор. После выполнения необходимых вычислений отображается следующее окно со схемой планшета:



В каждой лунке отображается соответствующее вычисленное значение. Вид окна со схемой планшета зависит от выбора элементов на панели управления, используемых для изменения параметров и настроек. Если требуется изменить метод, нажмите на вкладке **Редактировать метод**.



Если нажать на лунке правой кнопкой мыши, отображается следующее контекстное меню:

SM1	_10 SM1_14 SM1_18 SM1_2		
2. 0.1	Сводка Подробности Водаткировости		
SM1 1. 0.0	Изменить кинетически настройки Копировать кинетические настройки Вставить кинетические настройки		
SM1 2. 0	График: несколько меток График: кинетический анализ График: несколько меток кинетики График: ВЖФ кривых		
SM1	График: спектров График: серии разведений		
0.0	Выбор маскировать/показать Показать/Скрыть разметку		
SM1	12 SM1 16 SM1 20 SM1 2		

При выборе пункта Сводка отображается следующее окно с подробной информацией об определении и настройках выбранной лунки:

Лунка: С5 🛛 🔀		
	-	
😑 🛲 Схема метода		
🕀 🛲 Идентиф.		
🕀 📠 Псевдон. жидк		
Повторение		
<ul> <li>— Дазведение</li> </ul>		
🛨 🛲 Обработка данных для нескольких меток		
🕂 🚐 Формула(лы) преобразования		
— 🚐 Формула(лы) кинетического преобразования		
— 📠 Формула(лы) преобразования концентрации		
🗄 📠 Имена метода для стрипов		
🗄 🌐 Данные прибора		
🗄 🔆 Уменьшенные данные		
🕀 🔟 Средние значения		
🛨 🔏 Преобразованные данные		
🛨 🥫 Концентрации		
🗄 🔟 Статистика		
Влево Вверх Вправо Вниз		
Раскрыть все ОК	]	

Нажмите в окне со схемой планшета кнопку **Далее** для открытия диалогового окна **Сохранить как**, в котором можно ввести имя файла и комментарии. Для сохранения файла нажмите небольшую кнопку **Сохранить** в левой части окна. При этом можно продолжить работу с методом или рабочим пространством. Для сохранения файла и завершения работы мастера нажмите кнопку **Готово** в правом нижнем углу экрана. Снова открывается список мастеров.



### 6.2.6 Определение количественного анализа ELISA

#### 1. Вычитание значения бланка

#### Определения в Magellan

Выберите в панели управления элемент **Добавить новое преобразование**. Отображается диалоговое окно с запросом о необходимости определения **учета нулевого уровня**. После нажатия кнопки **Да** формула **Учет нулевого уровня** автоматически применяется в отношении всех лунок.

#### 2. Определение концентраций

## Определения в Magellan (Панель управления – Схема метода/ Конц.-, Разв.-, Эталон.-значения)

Выбранный идентификатор: ST Единица: УЕ/мл

ST1_1	5	(ST1_1Стандарт (калибратор) 1 для экспериментальной группы 1)
ST1_2	10	(ST1_2Стандарт (калибратор) 2 для экспериментальной группы 1)
ST1_3	20	(ST1_3Стандарт (калибратор) 3 для экспериментальной группы 1)
ST1_4	40	(ST1_4Стандарт (калибратор) 4 для экспериментальной группы 1)
ST1_5	80	(ST1_5Стандарт (калибратор) 5 для экспериментальной группы 1)
ST1_6	160	(ST1_6Стандарт (калибратор) 6 для экспериментальной группы 1)

#### 3. Определение калибровочной кривой

## Определения в Magellan (Панель управления – Концентрации/ Калибровочная кривая)

Входные данные:	учет нулевого уровня
Тип анализа	линейная регрессия
Ось Х	линейная
Ось Ү	линейная

#### 4. Определение граничных значений

## Определения в Magellan (Панель управления – Оценка данных/ Определение граничных значений)

Входные данные: средняя концентрация (УЕ/мл) Пределы: 22 18 положительный ≥ 22 > промежуточный ≥ 18 > отрицательный

Неконкурентный тест



#### 5. Контроль качества

## Определения в Magellan (Панель управления – Оценка данных/ Контроль качества)

Входные данные: единичная концентрация (УЕ/мл) Условия проверки 1 NC1\_1<8 Условия проверки 2 NC1\_2<8

NC1\_1....отрицательный контроль первого повторения для экспериментальной группы 1

NC1\_2.....отрицательный контроль второго повторения для экспериментальной группы 1



## 7. Чистка, техническое обслуживание и утилизация

## 7.1 Введение

В этой главе описаны следующие процедуры:

- чистка прибора;
- дезинфекция прибора;
- техническое обслуживание прибора;
- установка или замена фильтров в диске фильтров;
- утилизация.



#### ПРЕДУПРЕЖДЕНИЕ ПЕРЕД ВЫПОЛНЕНИЕМ ЛЮБЫХ РАБОТ ПО ЧИСТКЕ ИЛИ ТЕХНИЧЕСКОМУ ОБСЛУЖИВАНИЮ ПРИБОРА НЕОБХОДИМО СНЯТЬ МИКРОПЛАНШЕТ.



#### ПРЕДУПРЕЖДЕНИЕ

ПЕРЕД ВЫПОЛНЕНИЕМ РАБОТ ПО ЧИСТКЕ ИЛИ ДЕЗИНФЕКЦИИ ПРИБОРА ОТСОЕДИНИТЕ ЕГО ОТ ВНЕШНЕГО ИСТОЧНИКА ПИТАНИЯ.



#### ПРЕДОСТЕРЕЖЕНИЕ ЗАПРЕЩАЕТСЯ ВРУЧНУЮ ПЕРЕМЕЩАТЬ ТРАНСПОРТ ПЛАНШЕТА, ЕСЛИ ПРИБОР НЕ ВЫКЛЮЧЕН.

## 7.2 Чистка прибора



### ПРЕДУПРЕЖДЕНИЕ

ЧИСТКА ДОЛЖНА ПРОИЗВОДИТЬСЯ В ХОРОШО ПРОВЕТРИВАЕМОМ ПОМЕЩЕНИИ УПОЛНОМОЧЕННЫМ ПОДГОТОВЛЕННЫМ СПЕЦИАЛИСТОМ, ИСПОЛЬЗУЮЩИМ ОДНОРАЗОВЫЕ ПЕРЧАТКИ, ЗАЩИТНЫЕ ОЧКИ И ЗАЩИТНУЮ ОДЕЖДУ.

Очищайте корпус прибора и транспорт планшета только сухой или влажной тканью. При сильном загрязнении очистите прибор тканью, смоченной этиловым спиртом в концентрации не более 70% или раствором мягкого моющего средства (Microcide SQ или Decon 90). Протрите насухо бумажной безворсовой салфеткой.

Если на прибор попала какая-либо жидкость, ее следует немедленно удалить, чтобы предотвратить попадание жидкости в оптическую систему, которое может вызвать нарушение работоспособности прибора или сообщение об ошибке.



## 7.3 Дезинфекция прибора



#### ПРЕДУПРЕЖДЕНИЕ

ЕСЛИ ЖИДКОСТЬ, ПРОТЕКШАЯ НА ТРАНСПОРТ ПЛАНШЕТА, ПРЕДСТАВЛЯЕТ ИНФЕКЦИОННУЮ ОПАСНОСТЬ, СЛЕДУЕТ ПРОВЕСТИ ДЕЗИНФЕКЦИЮ СОГЛАСНО СООТВЕТСТВУЮЩИМ ГОСУДАРСТВЕННЫМ ЗАКОНАМ И НОРМАТИВАМ.

Все детали прибора, которые контактировали с биологическими образцами, образцами от пациентов, положительными контролями или биологически опасным материалом, должны рассматриваться как потенциально инфицированные участки.



#### ПРЕДУПРЕЖДЕНИЕ ПРОЦЕДУРА ДЕЗИНФЕКЦИИ И ДЕЗИНФИЦИРУЮЩИЕ СРЕДСТВА ДОЛЖНЫ СООТВЕТСТВОВАТЬ ТРЕБОВАНИЯМ ПРИМЕНИМЫХ ГОСУДАРСТВЕННЫХ ЗАКОНОВ И НОРМАТИВНЫХ АКТОВ.



#### ПРЕДУПРЕЖДЕНИЕ ОЧЕНЬ ВАЖНО ВЫПОЛНЯТЬ ТЩАТЕЛЬНУЮ ДЕЗИНФЕКЦИЮ ПРИБОРА ПЕРЕД ЕГО ВЫНОСОМ ИЗ ЛАБОРАТОРИИ ИЛИ ВЫПОЛНЕНИЕМ КАКОГО-ЛИБО ТЕХНИЧЕСКОГО ОБСЛУЖИВАНИЯ.

Перед возвращением прибора торговому представителю или в сервисный центр все внешние поверхности и транспорт планшета должны быть продезинфицированы. Организация, эксплуатирующая прибор, должна заполнить сертификат безопасности. При отсутствии сертификата безопасности прибор может быть не принят торговым представителем или сервисным центром или его могут задержать таможенные органы.

## 7.3.1 Растворы для дезинфекции

Дезинфекцию внешних поверхностей и транспорта прибора следует выполнять при помощи следующих растворов для дезинфекции поверхностей:

- Microcide SQ
- Decon 90
- 70 %-й раствор этилового спирта



ОПАСНОСТЬ ПОЖАРА И ВЗРЫВА! СПИРТЫ, ТАКИЕ КАК ЭТИЛОВЫЙ ИЛИ ИЗОПРОПИЛОВЫЙ, ЯВЛЯЮТСЯ ЛЕГКОВОСПЛАМЕНИМЫМИ И ПРИ НЕПРАВИЛЬНОМ ОБРАЩЕНИИ МОГУТ ВЫЗВАТЬ ВЗРЫВ ИЛИ ПОЖАР. ПРИ ВЫПОЛНЕНИИ РАБОТ С НИМИ СОБЛЮДАЙТЕ НАДЛЕЖАЩИЕ ПРАВИЛА ТЕХНИКИ БЕЗОПАСНОСТИ ДЛЯ ЛАБОРАТОРИЙ.

ПРЕДУПРЕЖДЕНИЕ





#### ПРЕДОСТЕРЕЖЕНИЕ ЗАПРЕЩАЕТСЯ ПРИМЕНЕНИЕ АЦЕТОНА, ПОСКОЛЬКУ ЭТО ПРИВОДИТ К ПОВРЕЖДЕНИЮ КРЫШЕК.

## 7.3.2 Процедура дезинфекции

Если лаборатория не использует специальную процедуру дезинфекции, для дезинфекции наружных поверхностей прибора и транспорта планшета следует применять описанную ниже процедуру.







## ПРЕДУПРЕЖДЕНИЕ

ДЕЗИНФЕКЦИЯ ДОЛЖНА ПРОИЗВОДИТЬСЯ В ХОРОШО ПРОВЕТРИВАЕМОМ ПОМЕЩЕНИИ УПОЛНОМОЧЕННЫМ ПОДГОТОВЛЕННЫМ СПЕЦИАЛИСТОМ, ИСПОЛЬЗУЮЩИМ ОДНОРАЗОВЫЕ ПЕРЧАТКИ, ЗАЩИТНЫЕ ОЧКИ И ЗАЩИТНУЮ ОДЕЖДУ.

#### ПРЕДОСТЕРЕЖЕНИЕ

СРЕДСТВО ДЕЗИНФЕКЦИИ ПОВЕРХНОСТИ МОЖЕТ ОТРИЦАТЕЛЬНО ПОВЛИЯТЬ НА РАБОТУ ПРИБОРА В СЛУЧАЕ ЕГО ПРИМЕНЕНИЯ ВНУТРИ ПРИБОРА ИЛИ СЛУЧАЙНОГО ПРОНИКНОВЕНИЯ ТУДА.

#### ПРЕДУПРЕЖДЕНИЕ

ВО ИЗБЕЖАНИЕ ОПАСНОСТИ ПОЖАРА ИЛИ ВЗРЫВА ПЕРЕД ДЕЗИНФЕКЦИЕЙ ПРИБОР СЛЕДУЕТ ОТСОЕДИНИТЬ ОТ ЭЛЕКТРОСЕТИ.

- 1. Наденьте защитные перчатки, защитные очки и защитную одежду.
- 2. Подготовьте подходящий контейнер для всех одноразовых материалов, используемых во время процедуры дезинфекции.
- 3. Разорвите связь между прибором и программным обеспечением и выключите прибор.



#### 7. Чистка, техническое обслуживание и утилизация

4. Осторожно выдвиньте транспорт планшета из прибора.



- Аккуратно нанесите раствор для дезинфекции на транспорт планшета согласно инструкциям производителя.
   Во избежание проникновения раствора внутрь прибора или загрязнения линз при задвижении транспорта планшета не наносите слишком много раствора.
- 6. По прошествии необходимого времени воздействия (согласно инструкциям производителя по применению) протрите транспорт планшета мягким бумажным полотенцем, смоченным мягким моющим средством или дистиллированной водой, чтобы удалить все остатки дезинфицирующего средства.
- 7. Осторожно задвиньте транспорт планшета в прибор.



- 8. Аккуратно нанесите раствор для дезинфекции на основание транспорта планшета.
- По прошествии необходимого времени воздействия протрите основание транспорта планшета мягким бумажным полотенцем, смоченным мягким моющим средством или дистиллированной водой, чтобы удалить все остатки дезинфицирующего средства.



- 10. Аккуратно нанесите раствор для дезинфекции на все наружные поверхности прибора.
- 11.По прошествии необходимого времени воздействия протрите прибор мягким бумажным полотенцем, смоченным мягким моющим средством или дистиллированной водой, чтобы удалить все остатки дезинфицирующего средства.
- 12.Протрите насухо наружные поверхности прибора мягким бумажным полотенцем.
- 13.Повторите процедуру дезинфекции для всех принадлежностей, перенесенных в другое место или возвращенных к прибору.
- 14.Уничтожьте контейнер с одноразовыми материалами в соответствии с применимыми государственными законами и нормативами.
- 15. Продезинфицируйте руки и вымойте их мягким моющим средством.

Перед возвратом прибора торговому представителю или в сервисный центр выполните следующее:

16. Упакуйте прибор и его принадлежности.

17.Составьте сертификат безопасности (см. ниже) и прикрепите его на упаковку в заметном месте.

### 7.3.3 Сертификат безопасности

Чтобы обеспечить безопасность и здоровье персонала сервисного центра, перед отправкой туда прибора для выполнения технического обслуживания или ремонта заполните бланк **Сертификата безопасности** (входящий в комплект поставки прибора) и прикрепите одну копию на верхней стороне упаковки возвращаемого прибора (в заметном месте!). Другую копию включите в сопроводительные документы.

Перед отправкой прибора эксплуатирующая организация обязана продезинфицировать его (см. 7.3.2 Процедура дезинфекции).

Дезинфекция должна производиться в хорошо проветриваемом помещении уполномоченным подготовленным специалистом, использующим неопудренные перчатки, защитные очки и защитную одежду.

Процедура дезинфекции должна производиться в соответствии с применимыми государственными, региональными и местными предписаниями.

При отсутствии сертификата безопасности прибор может быть не принят сервисным центром.

При необходимости можно запросить новую копию сертификата безопасности в сервисном центре Tecan.

# 7.4 План профилактического технического обслуживания фотометра INFINITE F50 PLUS

Рекомендуется выполнять следующие процедуры профилактического технического обслуживания.

**TECAN** 

### 7.4.1 Ежемесячно

Чистка корпуса и транспорта планшета с применением мягкого моющего средства (при необходимости чаще).



#### ПРЕДОСТЕРЕЖЕНИЕ ЗАПРЕЩАЕТСЯ ПРИМЕНЕНИЕ АЦЕТОНА, ПОСКОЛЬКУ ЭТО ПРИВОДИТ К ПОВРЕЖДЕНИЮ КРЫШЕК.

## 7.4.2 Каждые четыре года

Замена фильтров.

## 7.5 Замена и установка фильтров

Для выполнения установки или замены фильтров фотометр INFINITE F50 PLUS необходимо подсоединить к Magellan. В случае случайного нарушения в ходе этой процедуры связи компьютера и прибора необходимо выйти из Magellan и выключить прибор. Для возобновления процедуры выполните приведенные ниже действия. По их завершении восстановите соединение, для чего перезапустите устройство и Magellan и определите вновь установленные фильтры.



#### ПРЕДОСТЕРЕЖЕНИЕ

ПРИ ОБРАЩЕНИИ С ФИЛЬТРАМИ, СЛЕДИТЕ ЗА ТЕМ, ЧТОБЫ НЕ ПОЦАРАПАТЬ И НЕ ЗАГРЯЗНИТЬ ИХ ОТПЕЧАТКАМИ ПАЛЬЦЕВ ИЛИ ПЫЛЬЮ.



## 7.5.1 Замена фильтров

Для установки новых фильтров в стандартный диск фильтров или замены существующих фильтров выполните следующие действия:

- 1. В окне «Список мастеров» выберите Дополнительно.
- 2. Щелкните Управление прибором.
- 3. Щелкните Определение кассет фильтров.
- 4. Для запуска процедуры щелкните Замена фильтров.
- 5. Снимите с транспорта все микропланшеты!
- 6. Осторожно опрокиньте прибор назад и положите его на заднюю сторону дном к себе.
- Выверните оранжевые винты и снимите нижнюю крышку прибора.



8. Осторожно извлеките диск фильтров. Имейте в виду, что диск имеет магнитное крепление.



9. Поместите диск на чистую плоскую поверхность.



#### 7. Чистка, техническое обслуживание и утилизация

 Для выполнения замены фильтра извлеките фильтр из гнезда, пользуясь специальным монтажным инструментом.
 Для заказа этого инструмента и фильтров обратитесь к местному



- 11.Совместите монтажный инструмент с вырезом стопорного кольца. Проверните инструмент и извлеките стопорное кольцо из гнезда фильтра.
- 12. Проверните диск фильтров так, чтобы фильтр мог выскользнуть из гнезда. Чтобы не поцарапать фильтр, не выталкивайте его монтажным инструментом.
- 13. Установите в гнездо новый фильтр с учетом его ориентации. Действуйте осторожно, чтобы не поцарапать фильтр и не загрязнить его отпечатками пальцев.





Примечание Убедитесь в правильности установки фильтра.


#### 7. Чистка, техническое обслуживание и утилизация

14.Поместите стопорное кольцо на торец монтажного инструмента и поверните так, чтобы оно не соскользнуло.



- 15. Используя монтажный инструмент, вдавите стопорное кольцо в гнездо фильтра до достижения плотной посадки.
- 16. Проверните инструмент до совмещения его торца с разрывом стопорного кольца, после чего снимите инструмент.



- 17.Поместите диск фильтров обратно в гнездо диска и вдавите глубже, пока он не притянется магнитом.
- 18. Установите обратно крышку на дно прибора и притяните ее тремя оранжевыми винтами.
- 19.Поставьте прибор в нормальное положение.
- 20.Для завершения процедуры и запуска инициализации диска фильтров нажмите кнопку **ОК**.
- 21.Определите вновь установленный фильтр (процедура определения представлена в следующей главе).



### 7.5.2 Определение фильтров

В диалоговом окне Определение фильтра задайте новые длины волн в позициях, соответствующих замененным фильтрам.

Если фильтр установлен в новую позицию, активируйте эту позицию путем выбора ее флажка и введите соответствующую длину волны.

Для сохранения определений фильтров и запуска их инициализации нажмите кнопку **Сохранить**.

По завершении инициализации фильтров прибор готов к выполнению измерений.



Примечание Будьте внимательны! Перепутывание позиций и длин волн фильтров приведет к неверным результатам измерения.

# 7.6 Утилизация

### 7.6.1 Введение

Следуйте лабораторным процедурам по утилизации биологически опасных отходов в соответствии с государственными и местными предписаниями.

В этой главе приводятся инструкции по законной утилизации отходов, образующихся при использовании фотометра INFINITE F50 PLUS.



### ПРЕДОСТЕРЕЖЕНИЕ

СОБЛЮДАЙТЕ ВСЕ ГОСУДАРСТВЕННЫЕ, РЕГИОНАЛЬНЫЕ И МЕСТНЫЕ ПРЕДПИСАНИЯ ПО ОХРАНЕ ОКРУЖАЮЩЕЙ СРЕДЫ.

### 7.6.2 Утилизация упаковочных материалов

Упаковка прибора выполнена из материалов, пригодных для повторного использования. Если вы не намереваетесь использовать упаковку в будущем, например, для целей транспортировки или хранения, утилизируйте ее в соответствии с местными предписаниями.

## 7.6.3 Утилизация рабочих материалов

### ПРЕДУПРЕЖДЕНИЕ

В ПРОЦЕССЕ ЭКСПЛУАТАЦИИ ФОТОМЕТРА INFINITE F50 PLUS ОБРАЗУЮТСЯ ОТХОДЫ (МИКРОПЛАНШЕТ), ПРЕДСТАВЛЯЮЩИЕ БИОЛОГИЧЕСКУЮ ОПАСНОСТЬ.



ОБРАЩАЙТЕСЬ С МИКРОПЛАНШЕТАМИ, ДРУГИМИ ОДНОРАЗОВЫМИ МАТЕРИАЛАМИ И ВСЕМИ ИСПОЛЬЗУЕМЫМИ ВЕЩЕСТВАМИ В СООТВЕТСТВИИ С РЕКОМЕНДАЦИЯМИ ПО ЛАБОРАТОРНОЙ ПРАКТИКЕ.

ЗАПРОСИТЕ ИНФОРМАЦИЮ О СООТВЕТСТВУЮЩИХ ПРИЕМНЫХ ПУНКТАХ И УТВЕРЖДЕННЫХ МЕТОДАХ УТИЛИЗАЦИИ В ВАШЕЙ СТРАНЕ, РЕГИОНЕ ИЛИ ГОРОДЕ.



### 7.6.4 Утилизация прибора

В случае вопросов относительно утилизации прибора обратитесь в местный сервисный центр Tecan.

Степень загрязнения	2 (IEC/EN 61010-1)
Метод утилизации	Загрязненные отходы

X
1



(DIRECTIVE 2012/19/EU ON WASTE ELECTRICAL AND ELECTRONIC EQUIPMENT (WEEE)).

ВНИМАНИЕ ОТРИЦАТЕЛЬНОЕ ВЛИЯНИЕ НА ОКРУЖАЮЩУЮ СРЕДУ, СВЯЗАННОЕ С ДИРЕКТИВОЙ 2012/19/EU ПО ОБРАБОТКЕ ОТХОДОВ ЭЛЕКТРИЧЕСКОГО И ЭЛЕКТРОННОГО ОБОРУДОВАНИЯ

- НЕ ВЫБРАСЫВАЙТЕ ЭЛЕКТРИЧЕСКОЕ И ЭЛЕКТРОННОЕ ОБОРУДОВАНИЕ ВМЕСТЕ С НЕСОРТИРОВАННЫМ БЫТОВЫМ МУСОРОМ.
- СОБИРАЙТЕ ОТХОДЫ ЭЛЕКТРИЧЕСКОГО И ЭЛЕКТРОННОГО ОБОРУДОВАНИЯ ОТДЕЛЬНО.

### ПРЕДУПРЕЖДЕНИЕ

В ЗАВИСИМОСТИ ОТ ПРИМЕНЕНИЯ, НА ФОТОМЕТР INFINITE F50 PLUS МОГЛО ПОПАСТЬ ВЕЩЕСТВО, ПРЕДСТАВЛЯЮЩЕЕ БИОЛОГИЧЕСКУЮ ОПАСНОСТЬ.

- ОБЯЗАТЕЛЬНО ОБРАЩАЙТЕСЬ С НИМ В СООТВЕТСТВИИ С ПРИМЕНИМЫМИ СТАНДАРТАМИ И НОРМАТИВАМИ ПО БЕЗОПАСНОСТИ.
- ОБЯЗАТЕЛЬНО ДЕЗИНФИЦИРУЙТЕ ВСЕ ЧАСТИ ПЕРЕД УТИЛИЗАЦИЕЙ.

Для получения дополнительной информации о данном изделии обращайтесь в компанию Tecan:

Компания Tecan Austria GmbH

Untersbergstrasse 1A

A-5082 Grödig/Salzburg

AUSTRIA/EUROPE (ABCTPИЯ/EBPOПА)

Телефон: +43 6246 8933 444

Электронная почта: expertline-at@tecan.com.

www.tecan.com



# 8. Поиск неисправностей

## 8.1 Введение

Электронные функции прибора, а также измерения, операции и результаты, контролируются внутренним микропроцессором. При обнаружении микропроцессором неисправности или нештатного выполнения какой-либо процедуры на мониторе компьютера отображается сообщение об ошибке.

### 8.1.1 Таблица сообщений об ошибках и поиск неисправностей

В следующей таблице приведены краткие описания сообщений об ошибках и действий по поиску неисправностей.



Примечание При появлении сообщений об ошибках, не указанных в приведенной ниже таблице, обратитесь в местный сервисный центр Tecan.

Сообщение об ошибке	Описание	Поиск неисправности
Ошибка системы		
Lid Open Error (Ошибка! Открыта крышка)	В момент начала измерения открыта крышка	Закройте крышку и начните измерение заново
MTP Init Error (Ошибка инициализации микропланшета)	Не удалось инициализировать транспорт микропланшета	Аппаратная неисправность: неисправность электроники, обрыв ремня или механическая блокировка транспорта микропланшета
MTP lost steps (Потеря шагов микропланшета) abs( <i>steploss</i> ) > <i>max_steploss</i>	Потеря шагов микропланшета в ходе измерения <i>Steploss</i> : количество потерянных шагов <i>max_steploss</i> : допустимое количество потерянных шагов	Аппаратная неисправность: неисправность электронных компонентов, заедание механизма
Filter lost steps (Потеря шагов диска фильтров) abs( <i>steploss</i> ) > <i>max_steploss</i>	Потеря шагов диска фильтров в ходе измерения. <i>Steploss</i> : количество потерянных шагов <i>max_steploss</i> : допустимое количество потерянных шагов	Аппаратная неисправность: неисправность электронных компонентов, заедание механизма
USB timeout (Тайм-аут USB)	Тайм-аут в связи по USB	Ошибка системы – сообщите в сервисный центр



### 8. Поиск неисправностей

Сообщение об ошибке	Описание	Поиск неисправности
Ошибка системы		
Lamp Low! (Неисправность лампы) минимальное значение: <i>minimum</i> , максимальное значение: <i>maximum</i> Это сообщение отображается в микропрограммах версии V1.11 или более ранних версий	Измеренная интенсивность света находится вне нормативного диапазона между значениями <i>minimum</i> и <i>maximum</i>	Аппаратная неисправность: неисправность электронных компонентов, обрыв оптического волокна
Начиная с версии V1.12, сообщение «Lamp Low» заменено на сообщение «Prepare REF check»	Измеренная интенсивность света ниже ожидаемого диапазона между значениями <i>minimum</i> и <i>maximum</i>	Аппаратная неисправность: неисправность электронных компонентов, обрыв оптического волокна
Wavelength Not Available ! (Данная длина волны недоступна) Длина волны: <i>wavelength</i> нм	Фильтр с длиной волны <i>wavelength</i> не найден в диске фильтров	Ошибка системы – сообщите в сервисный центр
Channel Low! (Неисправность канала) Канал: <i>channel_nr</i> , минимальное значение: <i>minimum</i> , максимальное значение: <i>maximum</i>	Сигнал в канале <i>channel_nr</i> находится вне нормативного диапазона между значениями <i>minimum</i> и <i>maximum</i>	Аппаратная неисправность: неисправность электронных компонентов, обрыв оптического волокна
Неверная длина волны! Длина волны: <i>wavelength</i> нм	Длина волны фильтра находится вне диапазона длин волн белого и синего светодиодов	Ошибка системы – сообщите в сервисный центр
Lamp Overflow! (Переполнение по интенсивности света) Минимальное значение: <i>minimum</i> , максимальное значение: <i>maximum</i>	Сигнал в АЦП находится вне нормативного диапазона между значениями <i>minimum</i> и <i>maximum</i>	Аппаратная неисправность: неисправность электронных компонентов
Value Not Set (Значение не задано): <i>value -1</i>	Значение <i>value</i> не задано	Ошибка системы – сообщите в сервисный центр
Filter Init Error (Ошибка инициализации фильтра)	Не удалось инициализировать механизм перемещения диска фильтров	Аппаратная неисправность: неисправность электронных компонентов, механическая блокировка механизма перемещения диска фильтров

## 8.1.2 Переполнение

Если результат измерения оптической плотности не соответствует характеристикам прибора (> 4,0 OD), происходит переполнение и измеренное значение OD в данной лунке заменяется словом «Переполнение». Это действие осуществляется не прибором, а управляющим программным обеспечением.



### 8.1.3 Сбой электропитания

В этом разделе описаны последствия сбоя электропитания.

Сбой электропитания прибора, но не управляющего компьютера (например, когда компьютер запитан через ИБП): возможна потеря USB-связи между прибором и компьютером. Управляющее программное обеспечение создает сообщение об ошибке.

Одновременный сбой электропитания прибора и управляющего компьютера: необходимо перезагрузить компьютер. Результаты измерения недоступны.



# 9. Сокращения, товарные знаки и символы

# 9.1 Сокращения

В данном Руководстве по эксплуатации могут встретиться следующие сокращения, приводимые здесь в справочных целях.

А	ампер
AC	Alternating Current (переменный ток)
АЦП	аналого-цифровой преобразователь
ANSI/SBS	American National Standards Institute (Национальный институт стандартов CША)/Society for Biomolecular Screening (Общество биомолекулярного тестирования)
ASCII	American Standard Code for Information Interchange (американский стандартный код для обмена информацией)
ASTM	American Society for Testing and Material (Американское общество специалистов по испытаниям и материалам)
°C	градус Цельсия
CE	Маркировка о соответствии требованиям СЕ
CFR	Code of Federal Regulations (свод федеральных правил США)
СМ	сантиметр
DC	Direct Current (постоянный ток)
EC	European Community (Европейское сообщество)
ELISA	Enzyme-linked Immunosorbent Assay (твердофазный иммуноферментный анализ)
EN	European Norm (европейский стандарт)
°F	градус Фаренгейта
FDA	Food and Drug Administration (Управление США по контролю за продуктами и лекарствами)
Гц	герц
IEC	International Electrotechnical Commission (Международная электротехническая комиссия (МЭК))
ID	Identification (идентификация)
IFU	Instructions for Use (руководство по эксплуатации)
IQ	Installation Qualification (проверка установки)
IVD	In Vitro Diagnostics (диагностика «in vitro»)
IVDR	In Vitro Diagnostic Regulation (Регламент о медицинских изделиях для диагностики «in vitro») (EU) 2017/746
кг	килограмм
Л	литр
LED	Light Emitting Diode (светодиод)
ЛИС	лабораторная информационная система
МГ	миллиграмм
мл	миллилитр



#### 9. Сокращения, товарные знаки и символы

ММ	миллиметр
MTP	Microplate (микропланшет)
мкл	микролитр
NFM	Nonferrous Metal (цветной металл)
NIST	National Institute of Standards and Technology (Национальный институт стандартов и технологий США)
НМ	нанометр
NRTL	Nationally Recognized Testing Laboratory (национальная испытательная лаборатория)
OD	Optical Density (оптическая плотность (OD))
OQ	Operational Qualification (эксплуатационная квалификация)
PCB	Printed Circuit Board (Доска РСВ Печатная)
RF	Radio Frequency (радиочастота)
RoHS	Restriction of the Use of Certain Hazardous Substances (ограничение использования определенных вредных веществ)
SOP	Standard Operating Procedure (стандартная рабочая процедура)
USB	Universal Serial Bus (универсальная последовательная шина)
УЕ	Условные единицы (измерения)
ΤÜV	Technischer Überwachungsverein (Союз работников технического надзора (ФРГ))
В	вольт
BA	вольт-ампер
WEEE	Waste Electrical and Electronic Equipment (отходы электрического и электронного оборудования)

# 9.2 Торговые знаки

Следующие названия продуктов и любые зарегистрированные и незарегистрированные товарные знаки, упомянутые в настоящем документе, используются только для целей идентификации и остаются в исключительной собственности соответствующих владельцев:

- Magellan<sup>™</sup>, Infinite<sup>®</sup>, MultiCheck<sup>™</sup>, Tecan<sup>®</sup> и Tecan Logo являются торговыми знаками компании Tecan Group Ltd., Männedorf, Switzerland (Швейцария);
- Windows<sup>®</sup> и Excel<sup>®</sup> являются зарегистрированными товарными знаками корпорации Microsoft Corporation, Redmond, WA, USA (США);
- Pentium<sup>®</sup> и Atom<sup>™</sup> являются товарными знаками корпорации Intel Corporation, Santa Clara, CA, USA (США);
- Adobe<sup>®</sup> Reader<sup>®</sup> является зарегистрированным товарным знаком корпорации Adobe Systems Incorporated, Seattle, WA, USA (США)
- Microcide SQ<sup>™</sup> является товарным знаком корпорации Global Biotechnologies Inc., Portland, ME, USA (США);
- Decon 90<sup>TM</sup> является товарным знаком компании Decon Laboratories Ltd., Hove, East Sussex, UK (Великобритания).



# 9.3 Символы

	Производитель
$\sim$	Дата изготовления
CE	Сертификат СЕ
UK CA	Маркировка United Kingdom Conformity Assessed указывает на соответствие изделия применимым нормам и правилам Великобритании.
ĺ	Перед работой с прибором прочитайте Руководство по эксплуатации
IVD	Медицинское устройство диагностики «In vitro»
UDI	Unique Device Identification Символ UDI на этикетке идентифицирует носитель данных.
REF	Номер заказа
REF SN	Номер заказа Серийный номер
REF SN ●←	Номер заказа Серийный номер Символ USB
REF SN •••••• X	Номер заказа Серийный номер Символ USB Символ соблюдения требований по отходам электрического и электронного оборудования (WEEE)
REF SN • XX 	Номер заказа Серийный номер Символ USB Символ соблюдения требований по отходам электрического и электронного оборудования (WEEE) Символ RoHS, Китай
REF SN $\leftarrow \leftarrow \leftarrow$ $\leftarrow \leftarrow$ $\leftarrow$ $\leftarrow \leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$ $\leftarrow$	Номер заказа Серийный номер Символ USB Символ соблюдения требований по отходам электрического и электронного оборудования (WEEE) Символ RoHS, Китай

На приборе имеются следующие символы.



# Указатель

### М

Magellan 17
-------------

## S

Software	
Magellan	
Magellan Tracker	

### В

Включение прибора		16
Встряхивание	. 23, 3	33

## Д

Дезинфекция	65, 66
Процедура	67
Растворы	66
Сертификат безопасности	69
Действия Панель управления	23
Диск фильтров	35
Дополнительно Панель управления.	25

## 3

Замена фильтров	. 70,	71
Запрос пользователя		25

### И

Измерения конечной точки	27
Измерения Панель управления	23
Измерения с несколькими метками	28
Инкубация	25
Информационная панель	27

## К

Кинетика Панель управления	24
Кинетические измерения	29
Кинетический цикл.	24
Комментарии	25
Контроль качества	39

## Л

Лаб. оборудование Панель управления	ı 22
-------------------------------------	------

### М

Мастер	
Запуск измерения	19
Оценка результатов	20
Прикрепление подписи	20
Создание/редактирование метода	20
Создание/редактирование списка	
идентификаторов образцов	20
Список мастеров	19
Микропланшеты	38
Обеспечение наилучших характеристи	к 32

## 0

Ожидать (Время)	25
Определение измерений	27
Определение фильтров	74
Оптимизация эксплуатационных	
характеристик	31
Оптическая плотность	23

### Π

Панель рабочего процесса	. 26
Панель управления	. 22
Действия	. 23
Дополнительно	. 25
Измерения	. 23
Кинетика	. 24
Лаб. оборудование	. 22
Переполнение	. 78
Планшет	. 22
Поиск и устранение неисправностей	. 77
Прибор	
Описание	. 33
Принадлежности	. 38
Расположение	. 31
Технические характеристики	. 36
Пример применения	. 43
Количественный анализ ELISA	. 43
Программное обеспечение	. 17
Пользовательский интерфейс	. 19
Проверка установки	. 18
Список мастеров	. 19
Установка	. 17
Программные элементы	
Встряхивание	. 23
Запрос пользователя	. 25
Инкубация	. 25
Кинетический цикл	. 24
Комментарии	. 25
Ожидать (Время)	. 25
Оптическая плотность	. 23
Планшет	. 22
Создание отступов и разблокирование	. 30
Условия кинетики	. 24
Часть планшета	. 22
Профиль пользователя	. 11
Процедура самопроверки	. 39

### Ρ

Распаковка	
Порядок распаковки 1	13
Распаковка и осмотр 1	13
Редактор параметров измерения 2	21
с	

9
;



### Указатель

Сертификат безопасности	69
Символы	83
Сокращения	81
Сообщения об ошибках	77
Список мастеров	19

### Т

Техника безопасности	7
Технические характеристики	
Техническое обслуживание	. 65, 70
План технического обслуживания	70
Требования к окружающей среде	14
Требования к питанию	14
Требования к системе	15
У	

Условия кинетики24
--------------------

Утилизация	65, 74
Прибор	75
Рабочий материал	74
Упаковочный материал	74
ч	

Часть планшета	. 22
Чистка	. 65

### Э

Эксплуатационная квалификация	39
Проверка сходимости	40
Тест MultiCheck	39
Тест линейности	41
Тестирование микропланшета	39



# Сервисные центры Тесап

Если у вас появятся какие-либо вопросы или потребуется техническая поддержка для вашего изделия Тесап, обратитесь в местный сервисный центр компании. Контактные данные можно найти на веб-сайте <a href="http://www.tecan.com/">http://www.tecan.com/</a>.

Чтобы получить максимальную поддержку, перед обращением в сервисный центр подготовьте следующие сведения (указанные на паспортной табличке изделия):

- название модели;
- серийный номер (SN);
- программное обеспечение с указанием версии (если применимо);
- описание проблемы и лицо для связи;
- дата и время возникновения проблемы;
- предпринятые вами шаги по устранению проблемы;
- ваши контактные данные (телефон, факс, адрес электронной почты и т. д.).

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.