

# thermoscientific

# Vanquish

# Fluoreszenz-Detektoren

VC-D50, VC-D51, VF-D50, VF-D51

# Betriebsanleitung

4820.7901-DE Version 2.0 •

Dezember 2019



Copyright © 2019 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

#### Übersetzung der Original-Betriebsanleitung

Die Hardware-Beschreibungen in dieser Anleitung beziehen sich auf die Gerätetypen VC-D50-A, VC-D51-A, VF-D50-A, VF-D51-A.

#### Warenzeichen

Acrobat, Adobe, und Adobe Reader sind Warenzeichen von Adobe Systems Incorporated. Microsoft und Windows sind Warenzeichen von Microsoft Corporation. Torx ist ein Warenzeichen von Acument Intellectual Properties, LLC. Alle anderen Warenzeichen sind Eigentum von Thermo Fisher Scientific und ihren Tochtergesellschaften.

#### Haftungsausschluss

Dieses Dokument liegt den Produkten von Thermo Fisher Scientific Inc. beim Kauf bei und ist beim Betrieb des Produkts zu beachten. Dieses Dokument ist urheberrechtlich geschützt; jedes teilweise oder vollständige Vervielfältigen dieses Dokuments ist ohne vorherige schriftliche Genehmigung durch Thermo Fisher Scientific Inc. untersagt.

Das vorliegende Handbuch wurde nach bestem Wissen und Gewissen erstellt. Der Inhalt dieses Handbuchs kann jederzeit ohne Ankündigung in späteren Versionen geändert werden.

Thermo Fisher Scientific Inc. erhebt keinen Anspruch auf die Vollständigkeit, Korrektheit und Fehlerfreiheit dieses Dokuments. Thermo Fisher Scientific Inc. übernimmt keine Haftung für Fehler, Versäumnisse, Schäden oder Verluste, die aus dem Gebrauch dieses Dokuments entstehen, selbst wenn die Informationen in diesem Dokument genau befolgt werden.

Dieses Dokument ist nicht Teil des Kaufvertrages zwischen Thermo Fisher Scientific Inc. und einem Kunden. Dieses Dokument regelt oder ändert keine Geschäftsbedingungen. Bei widersprüchlichen Informationen zwischen den beiden Dokumenten gelten die Geschäftsbedingungen.

#### Nur Druckversion der Anleitung

Gedruckt in Deutschland auf 100% chlorfrei gebleichtem, hochweißen Papier, das in einem umweltfreundlichen Verfahren hergestellt wird. Das führt zu einem Papierprofil mit null CO<sub>2</sub>-Emissionen.

#### Adresse des Herstellers

Dionex Softron GmbH, Part of Thermo Fisher Scientific, Dornierstrasse 4, D-82110 Germering

# Kontaktinformationen

So nehmen Sie Kontakt mit uns auf:

### Bestellinformationen

Fragen zu Bestellinformationen oder zum Vertrieb der HPLC-Produkte beantwortet Ihnen gerne Ihre lokale Thermo Fisher Scientific-Vertriebsorganisation. Weitere Kontaktdaten finden Sie unter Contact Us auf http://www.thermofisher.com.

### Technische Unterstützung

Wenn Sie technische Unterstützung für HPLC-Produkte benötigen, kontaktieren Sie Ihren Thermo Fisher Scientific-Kundendienst vor Ort. Weitere Kontaktdaten finden Sie unter Contact Us auf http:// www.thermofisher.com.

# Inhalt

1	Ver	wendung dieser Anleitung 11					
	1.1	Über o	Über diese Anleitung				
	1.2	Konventionen					
		1.2.1 Sicherheitshinweise					
		1.2.2	Besondere Hinweise und zusätzliche Informationen	13			
		1.2.3	Typografische Konventionen	14			
	1.3	Refere	enzdokumentation	15			
2	Sicl	herhei	it	17			
	2.1	Sicher	heitssymbole und Signalwörter	18			
		2.1.1	Sicherheitssymbole und Signalwörter in dieser Anleitung	18			
		2.1.2	Beachtung dieser Anleitung	18			
		2.1.3	Sicherheitssymbole am Gerät	19			
		2.1.4	Typenschild	19			
	2.2	Verwe	endungszweck	20			
	2.3	Sicher	heitsmaßnahmen	21			
		2.3.1	Allgemeine Sicherheitsinformationen	21			
		2.3.2	Qualifikation des Personals	22			
		2.3.3	Persönliche Schutzausrüstung	22			
		2.3.4	Sicherheitsmaßnahmen beim Umgang mit elektrischen Geräten	23			
		2.3.5 Allgemeine Restrisiken		24			
		2.3.6	Verhalten im Notfall	26			
	2.4	Inforn	nationen zu Lösungsmitteln und Additiven	27			
		2.4.1	Allgemeine Kompatibilität	27			
		2.4.2	Erlaubte pH-Bereiche	27			
		2.4.3	Erlaubte Konzentrationen	28			
		2.4.4	Weitere Informationen	28			
	2.5	Inforn	nationen zur Konformität	29			
3	Üb	erblic	k über das Gerät	31			
	3.1	Aussta	attung des Detektors	32			
	3.2	Funkti	ionsprinzip	33			
	3.3	Innen	ansicht	36			
	3.4	Messz	zelle	37			
	3.5	5 Lampe					

	3.6	Erkennen von Undichtigkeiten (Leakerkennung) 40			
	3.7	Betrieb			
4	Aus	spacken 4			
	4.1	Auspa	acken	44	
	4.2	Liefer	umfang	46	
5	Inst	tallati	on	47	
	5.1	Sicher	heitshinweise für die Installation	48	
	5.2	Install	lieren des Gerätes	49	
	5.3	Anfor	derungen an den Aufstellungsort	51	
	5.0	5.3.1	Hinweise zur Stromversorgung	51	
		5.3.2	Netzkabel	52	
		5.3.3	Kondensation	52	
	5.4	Öffne	n des Innenraums	53	
	5.5	Aufba	uen der Hardware	54	
		5.5.1	Systemaufbau	54	
		5.5.2	Verbinden des Gerätes	57	
		5.5.3	Einbauen der Messzelle	. 62	
	5.6	Anschließen der Flussverbindungen		64	
		5.6.1	Allgemeine Informationen und Hinweise	64	
		5.6.2	Kapillar- und Schlauchführung durch das System	65	
		5.6.3	Installieren der Einsätze für die Zwischenwand	67	
		5.6.4	Anschließen von Fittingen, Kapillaren und Schläuchen	68	
		5.6.5	Flussverbindungen an der Messzelle	71	
		5.6.6	Ableiten von Flüssigkeiten in den Abfall	77	
	5.7	Messe	en des Drucks in der Messzelle	78	
		5.7.1	Messen des Rückdrucks der Waste-Leitung	80	
		5.7.2	Messen des Rückdrucks des Vanquish-Systems (ohne Messzelle)	81	
		5.7.3	Bestimmen Sie den Rückdruck der Transferkapillare und des zusätzlich Moduls (ohne Messzelle)	ien 83	
		5.7.4	Bestimmen des Rückdrucks der Messzelle	85	
		5.7.5	Berechnen des Drucks am Einlass der Messzelle	86	
	5.8	Einsch	nalten des Gerätes	88	
	5.9	Einricl	hten des Gerätes in der Software	89	
6	Bet	rieb		91	
	6.1	Einfüh	nrung in dieses Kapitel	92	

	6.2	Sicherheitshinweise zum Betrieb			
	6.3	Bedienelemente			
		6.3.1	Tastatur		
		6.3.2	Statusanzeigen		
	6.4	Ein- ur	nd Ausschalten		
	6.5	Vorber	reiten des Gerätes für den Betrieb 98		
	6.6	Hinwe	ise zum Gebrauch von Messzellen 100		
	6.7	Betrie	osmodi des Detektors 101		
		6.7.1	Single-Channel-Modus 101		
		6.7.2	Multi-Channel-Modus (Mehrkanalbetrieb, nur VF-Detektoren) 101		
		6.7.3	Zero-Order-Modus 102		
		6.7.4	Single Spectrum Scan (Erfassung von Einzelspektren) 103		
		6.7.5	FL Field Acquisition 105		
	6.8	Wichti	ge Einstellungen für den Betrieb 106		
	6.9	Optimi	ieren der Geräteleistung 108		
		6.9.1	Allgemeine Hinweise		
		6.9.2	Übersicht der Parameter zur Optimierung 109		
		6.9.3	Optimale Anregungs- und Emissionswellenlänge bestimmen 110		
		6.9.4	Sensitivity (Detektorempfindlichkeit) 111		
		6.9.5	Filterrad (Nur VF-Detektoren) 115		
		6.9.6	PMT (nur wenn zweiter PMT installiert ist) 116		
		6.9.7	Datenaufnahmerate und Ansprechzeit (Single-Channel-Modus) 117		
		6.9.8	Messzellen-Temperatur 118		
		6.9.9	Lampen-Modus 119		
		6.9.10	Basislinienverhalten 120		
	6.10	Außerl	betriebnahme des Gerätes 121		
		6.10.1	Kurzzeitige Außerbetriebnahme (Betriebsunterbrechung) 121		
		6.10.2	Langfristige Außerbetriebnahme 122		
		6.10.3	Wiederaufnehmen des Gerätebetriebs nach langfristiger		
			Auserbetriebnahme 125		
7	Wa	rtung	und Service 127		
	7.1	Einfüh	rung in Wartung und Service 128		
	7.2	Sicher	neitshinweise zu Wartung und Service 129		
	7.3	Allgem	eine Regeln für Wartung und Service131		
	7.4	Wartu	ng und Wartungsintervalle		
		7.4.1	Wartungszeitplan 132		
		7.4.2	Reinigen oder Dekontaminieren des Gerätes 133		

		7.4.3	Predictive Performance	135
		7.4.4	Überwachen der Lampenbetriebsdauer	136
	7.5	Durch	führung einer Wellenlängen-Kalibrierung	137
	7.6	Durch	führung einer Wellenlängenvalidierung	139
	7.7	Messz	elle	141
		7.7.1	Hinweise zum Gebrauch von Messzellen	141
		7.7.2	Ausbauen der Messzelle	142
		7.7.3	Reinigen der Messzelle	144
		7.7.4	Einbauen der Messzelle	145
	7.8	Tausch	nen der Sicherungen	146
	7.9	Aktual	isieren der Gerätefirmware	148
	7.10	Tausch	nen der Türen	150
	7.11	Transp	oortieren und Versenden des Gerätes	152
		7.11.1	Vorbereiten des Gerätes für den Transport	152
		7.11.2	Transportieren des Geräts an einen anderen Standort	153
		7.11.3	Versenden des Gerätes	154
	7.12	Tausch	nen des Moduleinschubs	156
		7.12.1	Entfernen des Moduleinschubs	156
		7.12.2	Zurückschicken des Moduleinschubs	158
		7.12.3	Installieren des Moduleinschubs	159
		7.12.4	Anschließen des Moduleinschubs	161
8	Feh	lersuc	he	163
•	0.1		nin a lafarmatian an ann Eablanacha	104
	8.1	Aligen	ieine informationen zur Feniersuche	164
	8.2	Meldu	ngen	166
	8.3	Störun	ngen beim Betrieb	171
		8.3.1	Beheben von Undichtigkeiten	171
		8.3.2	Weitere mögliche Gerätestörungen beim Betrieb	172
9	Spe	zifikat	tionen	175
	9.1	Leistur	ngsspezifikationen	176
		9.1.1	Detektor	176
		9.1.2	Messzellen	178
	9.2	Techni	ische Spezifikationen	179
10	Zub	ehör,	Ersatzteile und Verbrauchsmaterialien	181
	10.1	Allgem	neine Informationen	182

	10.2	Zubeh	örkit	183
	10.3	Optior	ales Zubehör	184
	10.4	Ersatzt	eile und Verbrauchsmaterialien	185
11	Anh	nang		187
	11.1	Inform	ationen zur Konformität	188
		11.1.1	Konformitätserklärungen	188
		11.1.2	WEEE-Konformität	189
		11.1.3	Einhaltung der FCC-Richtlinien	189
		11.1.4	Versionsgeschichte der Anleitung	189
	11.2	UV-Cu	toff-Wellenlängen von Lösungsmitteln	190
	11.3	Digital	I/O	191
	Ind	ex		195

# 1 Verwendung dieser Anleitung

Dieses Kapitel gibt einen Überblick über diese Anleitung und die in dieser Anleitung verwendeten Konventionen und macht Angaben zu Referenzdokumenten, die zusätzlich zu dieser Anleitung verfügbar sind.

# 1.1 Über diese Anleitung

Diese Anleitung beschreibt die funktionalen Elemente und das Funktionsprinzip Ihres Vanquish™-Geräts und enthält Anweisungen für Installation, Aufbau, Inbetriebnahme, Außerbetriebnahme, Betrieb, Wartung und Fehlersuche.

Dieses Handbuch soll Ihnen den gezielten Zugriff auf die Abschnitte ermöglichen, die für Sie als Anwender interessant sind. Verschaffen Sie sich einen Überblick über Ihr Gerät, indem Sie die Anleitung gründlich durchlesen.

Diese Anleitung enthält auch Sicherheitsinformationen, Vorsichtsmaßnahmen und spezielle Hinweise, die korrekt eingehalten werden müssen, um Personenschäden, die Beschädigung des Geräts oder den Verlust von Daten zu vermeiden.

Beachten Sie folgende Hinweise:

- Die Gerätekonfiguration kann variieren; daher müssen nicht alle Beschreibungen zwangsläufig auch auf Ihr Gerät zutreffen.
- Bezieht sich eine Beschreibung nur auf ein Modell oder eine Variante, so wird das Modell oder die Variante namentlich genannt.
- Die Abbildungen in dieser Betriebsanleitung dienen dem grundlegenden Verständnis. Sie können vom eigentlichen Modell des Geräts oder der Komponente abweichen. Dies hat jedoch keinen Einfluss auf die Beschreibungen. Aus den Abbildungen in dieser Anleitung können keine Ansprüche hergeleitet werden.

Den Beschreibungen in dieser Anleitung liegt die Annahme zugrunde, dass das Gerät als Teil des Vanquish-Systemturms installiert wird. Ist dies nicht der Fall, wird zusätzliche Hardware benötigt; diese muss separat bestellt werden. Die Informationen in dieser Anleitung gelten entsprechend.

# 1.2 Konventionen

Dieser Abschnitt beschreibt die Konventionen, die für diese Anleitung gelten.

### 1.2.1 Sicherheitshinweise

Die Sicherheitshinweise oder Vorsichtsmaßnahmen in dieser Anleitung erscheinen wie folgt:

- Sicherheitshinweise oder Vorsichtsmaßnahmen, die für die gesamte Anleitung und alle in dieser Anleitung enthaltenen Anweisungen gelten, finden Sie im Kapitel Sicherheit.
- Sicherheitshinweise oder Vorsichtsmaßnahmen, die für einen ganzen Abschnitt oder mehrere in einem Abschnitt enthaltene Anweisungen gelten, finden Sie am Anfang des Abschnitts, für den sie gelten.
- Sicherheitshinweise, die nur f
  ür einen bestimmten Abschnitt oder eine bestimmte Anweisung gelten, befinden sich in dem jeweiligen Abschnitt oder in der Anweisung, f
  ür die sie gelten. Sie heben sich vom restlichen Text ab.

Sicherheitshinweise beginnen meist mit einem Gefahrensymbol und/ oder einem Signalwort. Das Signalwort erscheint in Großbuchstaben und fett gedruckt.

Stellen Sie sicher, dass Sie alle in dieser Anleitung enthaltenen Sicherheitshinweise verstehen und befolgen.

### **1.2.2** Besondere Hinweise und zusätzliche Informationen

Spezielle Hinweise und zusätzliche Informationen in dieser Anleitung heben sich vom restlichen Text ab. Sie erscheinen mit Rahmen und sind entsprechend bezeichnet. Die Bezeichnung erscheint in Großbuchstaben und fett gedruckt.

### ACHTUNG

Kennzeichnet Informationen, die Ihnen helfen, Schäden am Gerät oder ungültige Testergebnisse zu vermeiden.

**TIPP** Kennzeichnet Informationen von allgemeinem Interesse oder hilfreiche Informationen, die Ihnen eine Aufgabe erleichtern oder Ihnen helfen können, die Leistung des Geräts zu optimieren.

### 1.2.3 Typografische Konventionen

Für die Beschreibungen in dieser Anleitung gelten die folgenden typographischen Konventionen:

### Dateneingabe und Datenausgabe

- Folgende Texte erscheinen **fett** gedruckt:
  - Eingaben, die Sie über die Tastatur vornehmen oder mit der Maus auswählen
  - Schaltflächen, die Sie auf dem Bildschirm anklicken
  - Befehle, die Sie über die Tastatur eingeben
  - Bezeichnungen, zum Beispiel von Dialogfeldern, Properties und Parametern
- Zur besseren Übersichtlichkeit werden lange Formulierungen und Dateipfade in gekürzter Form verwendet, zum Beispiel: Klicken Sie Start > Alle Programme > Thermo Chromeleon 7 > Services Manager > Start Instrument Controller.

#### Referenzen und Meldungen

- Verweise auf zusätzliche Dokumente erscheinen kursiv.
- Meldungen, die auf dem Bildschirm angezeigt werden, erscheinen in Anführungszeichen.

#### Perspektive

Wenn nicht anders angegeben, beziehen sich die Ausdrücke *links* und *rechts* in dieser Anleitung immer auf die Perspektive einer Person, die direkt vor dem Gerät steht.

#### Besonders wichtige Begriffe

Besonders wichtige Begriffe im Text erscheinen kursiv.

#### Elektronische Version der Betriebsanleitung (PDF)

Die elektronische Version (PDF) der Anleitung enthält zahlreiche Verweise, auf die Sie klicken können, um innerhalb der Anleitung zu navigieren. Dazu gehören:

- Überschriften im Inhaltsverzeichnis
- Indexeinträge
- Querverweise (blaue Schrift), zum Beispiel auf Abschnitte und Abbildungen

# 1.3 Referenzdokumentation

Zusätzlich zu dieser Betriebsanleitung stehen weitere Referenzdokumente zur Verfügung.

#### Hardware-Dokumentation

Zu den Hardware-Dokumenten gehören unter anderem:

- Betriebsanleitungen zu den anderen Modulen des Vanquish-Systems Eine gedruckte Version der Anleitung liegt dem Gerät bei der Lieferung bei.
- Betriebsanleitung für das Vanquish-System Eine gedruckte Version der Anleitung liegt der Lieferung des Sockels und Solvent Racks des Vanquish-Systems bei.
- Bedienungsanleitung zur Instrument Installation Qualification

**TIPP** Diese Anleitungen sind elektronisch als PDF-Dateien (Portable Document Format) verfügbar. Um die PDF-Dateien öffnen und lesen zu können, werden Adobe™ Reader™ oder Adobe™ Acrobat™ benötigt.

### Software-Dokumentation

Folgende Software-Dokumentation ist verfügbar:

 Chromeleon<sup>™</sup>-Hilfe und Anwenderdokumentation
 Die Chromeleon-Hilfe bietet umfangreiche Informationen und ausführliches Referenzmaterial zu allen Aspekten der Software.

Zusätzlich steht folgende Dokumentation zur Verfügung (abhängig von der Software-Version):

- Installationsanleitung Grundlegende Informationen zur Geräteinstallation und Gerätekonfiguration finden Sie im Installation Guide.
- Instrument Configuration Manager-Hilfe Spezifische Informationen zu einzelnen Geräten finden Sie in der Instrument Configuration Manager-Hilfe. In Chromeleon 7 werden Geräte als 'Module' bezeichnet.
- Quick Start Guide
   Der Quick Start Guide beschreibt die wichtigsten Elemente der
   Benutzeroberfläche und führt Sie schrittweise durch die wichtigsten
   Arbeitsabläufe.

• *Reference Card* Die *Reference Card* beschreibt die wichtigsten Arbeitsabläufe in Kurzform.

**TIPP** Die *Chromeleon-Hilfe* und Anwenderdokumentation werden mit der Software ausgeliefert.

Weitere Dokumente

Beachten Sie auch die Anwenderdokumentation, die von den Herstellern der Drittanbieter-Komponenten und Drittanbieter-Substanzen zur Verfügung gestellt wird, zum Beispiel Sicherheitsdatenblätter (SDB).

# 2 Sicherheit

In diesem Kapitel finden Sie allgemeine und spezifische Sicherheitsinformationen sowie Informationen zum Verwendungszweck des Geräts.

# 2.1 Sicherheitssymbole und Signalwörter

### 2.1.1 Sicherheitssymbole und Signalwörter in dieser Anleitung

Diese Anleitung enthält Sicherheitshinweise zum Schutz von Personen, welche das Gerät betreiben.

Die folgenden Sicherheitssymbole und Signalwörter werden in dieser Anleitung verwendet:



Beachten Sie stets die Sicherheitsinformationen. Fahren Sie erst dann mit den Arbeiten fort, wenn Sie die Informationen vollständig verstanden und die Folgen Ihres Handelns bedacht haben.



### VORSICHT

Kennzeichnet eine gefährliche Situation, die, wenn sie nicht vermieden wird, zu kleinen oder leichten Verletzungen führen kann.



### WARNUNG

Kennzeichnet eine gefährliche Situation, die, wenn sie nicht vermieden wird, zu ernsthaften Verletzungen führen kann.

### 2.1.2 Beachtung dieser Anleitung

Beachten Sie Folgendes:

- Lesen Sie die Anleitung aufmerksam durch, bevor Sie das Gerät installieren oder betreiben, so dass Sie mit dem Gerät und der Anleitung vertraut sind. Die Anleitung enthält wichtige Informationen zur Sicherheit der Anwender sowie zu Gebrauch und Wartung des Geräts.
- Bewahren Sie diese Anleitung stets in der N\u00e4he des Ger\u00e4ts auf, damit sie bei Bedarf schnell zur Hand ist.
- Bewahren Sie diese Anleitung auf und geben Sie diese an nachfolgende Anwender weiter.



Lesen, verstehen und beachten Sie alle Sicherheitshinweise und Vorsichtsmaßnahmen in dieser Anleitung.

## 2.1.3 Sicherheitssymbole am Gerät

In der Tabelle sind die Sicherheitssymbole aufgeführt, die am Gerät oder auf Aufklebern am Gerät angebracht sind. Befolgen Sie die Sicherheitshinweise in dieser Anleitung, um eine Verletzungsgefahr für den Bediener und/oder Schäden am Gerät zu vermeiden.

Symbol	Beschreibung
	Weist auf eine mögliche Gefährdung hin. Beachten Sie die Informationen in dieser Anleitung, um eine Verletzungsgefahr von Personen und/oder Schäden am Gerät zu vermeiden.
0	Stromversorgung eingeschaltet Stromversorgung ausgeschaltet
~	Weist auf Wechselstrom hin.
	Weist darauf hin, dass sich die Oberfläche während des Betriebs erhitzt. Berühren Sie diese Oberflächen nicht, solange diese noch heiß sind.

### 2.1.4 Typenschild

Das Typenschild ist auf dem Gerät bei den elektrischen Anschlüssen angebracht. Das Typenschild gibt Auskunft über Seriennummer, Teilenummer, Modulnamen, Revisionsnummer (wenn vorhanden), Netzspannung und Nennleistung der Sicherungen und nennt die Herstelleradresse.

**TIPP** Ein weiteres Typenschild an der Leckage-Wanne des Geräts gibt Auskunft über den Modulnamen, die Seriennummer, die Teilenummer und die Revisionsnummer (wenn vorhanden). Für die Kommunikation mit Thermo Fisher Scientific benötigen Sie zur leichteren Identifizierung Ihres Gerätes die Informationen dieses Typenschilds.

# 2.2 Verwendungszweck

Das Gerät ist zur Verwendung im Vanquish-System gedacht.

Das Vanquish-System ist zur Analyse von Verbindungsgemischen in Probenlösungen gedacht.

Das Gerät darf nur von geschultem und qualifiziertem Personal in einer Laborumgebung betrieben werden.

Das Gerät und das Vanquish-System wurden ausschließlich als allgemeines Laborgerät (GLE = General Laboratory Equipment) entwickelt.

Sie sind nicht für den Einsatz in diagnostischen Verfahren gedacht.

#### Laborpraxis

Thermo Fisher Scientific empfiehlt, dass sich das Labor, welches das Vanquish-System betreibt, an die Richtlinien der Guten Laborpraxis für LC-Analysen hält. Dazu gehört unter anderem:

- Verwendung geeigneter Standards
- Regelmäßiges Kalibrieren
- Festlegung und Einhaltung von Grenzwerten f
  ür die Mindesthaltbarkeit aller mit dem System verwendeten Verbrauchsmaterialien
- Betrieb des Systems entsprechend der verifizierten und validierten laboreigenen Testprozedur

# 2.3 Sicherheitsmaßnahmen

## 2.3.1 Allgemeine Sicherheitsinformationen

Alle Anwender müssen zu jeder Zeit während Installation, Betrieb, Fehlerbehebung, Wartung, Außerbetriebnahme und Transport des Geräts die allgemeinen Sicherheitsinformationen in diesem Abschnitt sowie alle anderen in dieser Anleitung aufgeführten spezifischen Sicherheitshinweise und Vorsichtsmaßnahmen beachten.



Wenn das Gerät nicht entsprechend den Angaben von Thermo Fisher Scientific eingesetzt wird, kann der durch das Gerät bereitgestellte Schutz beeinträchtigt werden. Beachten Sie Folgendes:

- Betreiben Sie das Gerät nur innerhalb der technischen Spezifikationen.
- Verwenden Sie ausschließlich Ersatzteile und zusätzliche Komponenten, Optionen und Peripheriegeräte, die von Thermo Fisher Scientific ausdrücklich für das Gerät autorisiert und freigegeben sind.
- Führen Sie nur die Arbeiten durch, die in dieser Betriebsanleitung und in weiteren Dokumenten für das Gerät beschrieben sind. Folgen Sie allen Anweisungen Schritt für Schritt und verwenden Sie die in der Anleitung empfohlenen Werkzeuge.
- Öffnen Sie das Gehäuse des Geräts und anderer Komponenten nur, wenn Sie in dieser Anleitung dazu ausdrücklich aufgefordert werden.
- Für Schäden materieller oder immaterieller Art, die sich aus einer nicht bestimmungsgemäßen Verwendung oder unsachgemäßen Anwendung des Geräts ergeben, kann Thermo Fisher Scientific keine Haftung übernehmen. Fragen zur bestimmungsgemäßen Verwendung beantwortet Ihnen Thermo Fisher Scientific gern.

### Sicherheitsstandard

Das Gerät hat Schutzklasse I (mit einem Schutzleiter verbunden). Das Gerät wurde nach internationalen Sicherheitsstandards hergestellt und getestet.

## 2.3.2 Qualifikation des Personals

Beachten Sie die folgenden Informationen zu den Qualifikationen, die Personen besitzen müssen, welche das Gerät installieren und/oder bedienen.



Die Installation des Geräts und die Herstellung der elektrischen Verbindungen müssen durch geschultes Personal und entsprechend der geltenden Vorschriften erfolgen.

- Thermo Fisher Scientific empfiehlt, die Installation stets von Service-Personal durchführen zu lassen, das von Thermo Fisher Scientific entsprechend zertifiziert wurde (im Folgenden kurz als Thermo Fisher Scientific-Servicetechniker bezeichnet).
- Wenn Installation und Aufbau des Moduls durch eine andere Person als einen Thermo Fisher Scientific-Servicetechniker erfolgen, trägt diejenige Person die Verantwortung dafür, dass die Sicherheit von Modul und System gewährleistet ist.



### **Allgemeiner Betrieb**

Das Gerät darf nur von geschultem und qualifiziertem Personal in einer Laborumgebung betrieben werden.

Alle Anwender müssen die Gefahren kennen, die von dem Gerät und den verwendeten Substanzen ausgehen. Alle Anwender sollten die relevanten Sicherheitsdatenblätter (SDB) beachten.

### 2.3.3 Persönliche Schutzausrüstung

Tragen Sie persönliche Schutzausrüstung und folgen Sie der Guten Laborpraxis, um sich vor Gefahrstoffen zu schützen. Dabei hängt die passende Schutzausrüstung von der Gefahr ab. Informationen zu den Gefahren und der erforderlichen Schutzausrüstung der Substanzen, mit denen Sie umgehen, entnehmen Sie bitte dem Sicherheitsdatenblatt des Herstellers.



In der Nähe Ihres Arbeitsplatzes sollten sich eine Einrichtung zum Spülen der Augen und ein Spülbecken befinden. Falls die Substanz in Kontakt mit Ihren Augen oder Ihrer Haut kommt, waschen Sie die betroffenen Stellen mit Wasser ab und nehmen Sie sofort ärztliche Hilfe in Anspruch.

Schutzkleidung	
	Tragen Sie zum Schutz vor Chemikalienspritzern, gefährlichen Flüssigkeiten oder anderer Kontamination angemessene Schutzkleidung, zum Beispiel einen Laborkittel.
Augenschutz	
	Tragen Sie zum Schutz der Augen geeigneten Augenschutz, zum Beispiel eine Schutzbrille mit Seitenschutz. Besteht ein Risiko spritzender Flüssigkeiten, ist eine Vollsichtschutzbrille (Korbbrille) erforderlich.
Handschuhe	
	Tragen Sie zum Schutz vor gefährlichen Flüssigkeiten und zum Schutz vor Verletzungen während Wartungs- oder Servicearbeiten geeignete Schutzhandschuhe.

### 2.3.4 Sicherheitsmaßnahmen beim Umgang mit elektrischen Geräten



### WARNUNG—Stromschlag oder Schäden am Gerät

Im Gerät treten hohe Spannungen auf, die zu einem Stromschlag führen oder Schäden am Gerät verursachen können.

- Führen Sie keine Veränderungen an den elektrischen Anschlüssen oder Erdungsanschlüssen durch.
- Wenn Sie Schäden an der Elektrik vermuten, ziehen Sie den Netzstecker und wenden Sie sich an den Thermo Fisher Scientific-Kundendienst.
- Öffnen Sie nicht das Gehäuse oder entfernen Sie keine Schutzabdeckungen, es sei denn, Sie werden in dieser Anleitung dazu ausdrücklich aufgefordert.
- Stellen Sie keine Flüssigkeitsbehälter auf dem Gerät ab. Auslaufende Flüssigkeit könnte in das Gerät gelangen und in Kontakt mit elektronischen Komponenten kommen, und so einen Kurzschluss auslösen. Stellen Sie Flüssigkeitsbehälter stattdessen in das Solvent Rack des Vanquish-Systems.

### 2.3.5 Allgemeine Restrisiken

Beachten Sie die folgenden allgemeinen Restrisiken, wenn Sie mit dem Gerät arbeiten:



### WARNUNG—Gefährliche Substanzen

Lösungsmittel, mobile Phasen, Proben und Reagenzien können giftige, krebserregende, erbgutschädigende, infektiöse oder anderweitig schädliche Substanzen enthalten. Der Umgang mit diesen Substanzen kann Gesundheits- und Sicherheitsrisiken darstellen.

- Vergewissern Sie sich, dass Sie die Eigenschaften aller von Ihnen eingesetzten Substanzen kennen. Vermeiden Sie den Kontakt mit schädlichen Substanzen. Behandeln Sie Substanzen im Zweifelsfall wie eine gesundheitsschädliche Substanz.
- Tragen Sie persönliche Schutzausrüstung je nach Gefahr und folgen Sie der Guten Laborpraxis.
- Verwenden Sie nur die Substanzmengen, die mindestens für die Probenanalyse erforderlich sind.
- Betreiben Sie das Gerät nicht in einer brandgefährdeten Umgebung.
- Vermeiden Sie die Ansammlung schädlicher Substanzen. Stellen Sie sicher, dass der Aufstellungsort gut belüftet ist.
- Entsorgen Sie Abfälle gesundheitsschädlicher Substanzen umweltgerecht und entsprechend der lokalen Bestimmungen. Halten Sie bei der Entsorgung der Abfälle ein geregeltes und genehmigtes Verfahren ein.



### WARNUNG—Biogefährdung

Biologisch gefährliches Material, zum Beispiel Mikroorganismen, Zellkulturen, Gewebe, Körperflüssigkeiten und andere biologische Stoffe können ansteckende Krankheiten übertragen. So vermeiden Sie Infektionen durch biologische Stoffe:

- Behandeln Sie alle biologischen Substanzen als potentiell infektiös.
- Tragen Sie persönliche Schutzausrüstung je nach Gefahr und folgen Sie der Guten Laborpraxis.
- Entsorgen Sie Abfälle von Biogefahrstoffen umweltgerecht und entsprechend der lokalen Bestimmungen. Halten Sie bei der Entsorgung der Abfälle ein geregeltes und genehmigtes Verfahren ein.



#### WARNUNG—Selbstentzündung von Lösungsmitteln

Lösungsmittel, deren Selbstentzündungstemperatur unter 150 °C liegt, können sich beim Kontakt mit heißen Oberflächen (zum Beispiel, aufgrund von Undichtigkeiten im Chromatographie-System) selbst entzünden.

Vermeiden Sie die Verwendung derartiger Lösungsmittel.



#### WARNUNG—Gefährliche Dämpfe

Mobile Phasen und Proben können flüchtige oder brennbare Lösungsmittel enthalten. Der Umgang mit diesen Substanzen kann Gesundheits- und Sicherheitsrisiken darstellen.

- Vermeiden Sie die Ansammlung dieser Substanzen. Stellen Sie sicher, dass der Aufstellungsort gut belüftet ist.
- Vermeiden Sie offenes Feuer und Funken.
- Betreiben Sie das Gerät nicht in einer Umgebung mit brennbaren Gasen und Dämpfen.



#### VORSICHT—Austreten von Gefahrstoffen aus PEEK-Kapillaren

Im Vanquish-System können Kapillaren aus PEEK verwendet sein. Wenn PEEK-Kapillaren aufquellen oder von Säuren angegriffen werden, können sie undicht werden oder bersten. Bei Kontakt mit einigen Chemikalien, wie zum Beispiel Trichlormethan (CHCl<sub>3</sub>), Dimethylsulfoxid (DMSO) oder Tetrahydrofuran (THF) kann es zum Aufquellen des PEEKs kommen. Konzentrierte Säuren wie Schwefel- und Salpetersäure oder ein Gemisch aus Hexan, Ethylacetat und Methanol können PEEK angreifen.

- Das Aufquellen oder der Kontakt stellen bei kurzen Spülzyklen jedoch kein Problem dar.
- Weitere Informationen zur chemischen Beständigkeit von PEEK können Sie der technischen Literatur entnehmen.



#### VORSICHT—Allergische Reaktion

Einige Kapillaren im Vanquish-System sind aus der Nickel-Kobalt-Legierung MP35N™ gefertigt. Hautkontakt mit diesem Material kann bei Personen, die gegen Nickel/Kobalt empfindlich sind, gegebenenfalls eine allergische Reaktion hervorrufen.



### VORSICHT—Funkenbildung durch elektrostatische Entladung

Lösungsmittel, die durch Kapillaren fließen, können sich selbsttätig statisch aufladen. Dieser Effekt tritt insbesondere in isolierenden Kapillaren und bei nicht-leitenden Lösungsmitteln (beispielsweise reinem Acetonitril) auf. Elektrostatische Entladung kann zu Funkenbildung führen und eine Brandgefahr darstellen.

Vermeiden Sie die Entstehung von statischer Elektrizität im Bereich des Chromatographie-Systems.

#### 2.3.6 Verhalten im Notfall



### WARNUNG—Sicherheitsgefährdung

Trennen Sie im Notfall das Gerät vom Stromnetz.

# 2.4 Informationen zu Lösungsmitteln und Additiven

### 2.4.1 Allgemeine Kompatibilität

Beachten Sie im Hinblick auf eine optimale Funktionalität des Vanquish-Systems die folgenden Empfehlungen zur Verwendung von Lösungsmitteln und Additiven:

- Zusammen mit dem System dürfen ausschließlich Reversed-Phasekompatible (RP) Lösungsmittel und Additive verwendet werden.
- Verwenden Sie nur Lösungsmittel und Additive, die mit allen Teilen im Flussweg kompatibel sind.

**TIPP** In einem Vanquish Core-System dürfen Normal-Phasen-kompatible (NP) Lösungsmittel und Additive verwendet werden, wenn die VC-Pumpen und VC-Autosampler mit den Komponenten aus dem Normal-Phasen-Kit (NP-Kit) modifiziert sind. Informieren Sie sich in den *Betriebsanleitungen* zu Pumpe und Autosampler.

### 2.4.2 Erlaubte pH-Bereiche

Zulässige pH-Bereiche (Standard-Systemkonfiguration):

System (Standard- konfiguration)	Erlaubte pH- Bereiche	Bemerkungen		
Vanquish Core	1-13	• pH-Werte von 2 oder weniger: Die		
Vanquish Horizon	2-12	Anwendungsdauer sollte möglichst kurz sein. Spülen Sie das System nach entsprechenden Anwendungen gründlich.		
Vanquish Flex		• <i>pH-Werte über 9,5 mit optischen Detektoren</i> : Vermeiden Sie die Verwendung von mobilen Phasen mit einem pH-Wert über 9,5 mit optischen Detektoren. Dies kann die Funktionalität und optische Leistung der Messzelle im Detektor beeinträchtigen.		
		<ul> <li>pH-Werte über 12: Können die elektrochemische Detektion beeinträchtigen. Entfernen Sie daher den Detektor fluidisch aus dem System, bevor hochalkalische Lösungsmittel zum Spülen des Systems verwendet werden.</li> </ul>		
		• Ammoniumhydroxid-haltige mobile Phasen: In seltenen Fällen wurde bei Reversed-Phase- (UHMW-PE) Kolbendichtungen eine verkürzte Lebensdauer beobachtet, wenn sie über längere Zeit ammoniumhydroxid-haltigen mobilen Phasen mit hohem pH-Wert ausgesetzt waren.		

## 2.4.3 Erlaubte Konzentrationen

System (Standard- konfiguration)	Chlorid	Puffer	Bemerkungen
Vanquish Core	0,1 mol/L oder weniger	1 mol/L oder weniger	<ul> <li>Hohe Chlorid-Konzentration: Die Anwendungsdauer sollte möglichst kurz sein. Spülen Sie</li> </ul>
Vanquish Horizon Vanquish Flex	1 mol/L oder weniger	-	<ul> <li>das System nach entsprechenden Anwendungen gründlich.</li> <li>Ammoniumhydroxid-haltige mobile Phasen: In seltenen Fällen wurde bei Reversed- Phase- (UHMW-PE) Kolbendichtungen eine verkürzte Lebensdauer beobachtet, wenn sie über längere Zeit ammoniumhydroxid-haltigen mobilen Phasen mit hohem pH- Wert ausgesetzt waren.</li> </ul>

Erlaubte Konzentrationen (Standard-Systemkonfiguration):

### 2.4.4 Weitere Informationen

- Genauere Informationen zu den Materialien, die im Flussweg des Geräts verwendet sind, finden Sie im Kapitel *Spezifikationen* in dieser Anleitung. Informationen zu den Materialien, die im Flussweg der anderen Module im Vanquish-System verwendet werden, finden Sie im Kapitel *Spezifikationen* in der *Betriebsanleitung* der entsprechenden Module.
- Beachten Sie alle spezifischen Empfehlungen in anderen Abschnitten dieser Anleitung. Informieren Sie sich auch in den Betriebsanleitungen aller Module des Vanquish-Systems. Darin finden Sie gegebenenfalls weitere Richtlinien und Informationen.
- Beachten Sie die allgemeinen Richtlinien und Empfehlungen zur Verwendung von Lösungsmitteln und Additiven im Chromatographie-System. Informieren Sie sich dazu in Verwenden von Lösungsmitteln und Additiven in der Vanquish-System-Betriebsanleitung.

# 2.5 Informationen zur Konformität

Thermo Fisher Scientific führt umfassende Tests und Beurteilungen seiner Produkte durch, um die vollständige Einhaltung anwendbarer nationaler und internationaler Bestimmungen zu gewährleisten. Das Gerät erfüllt bei Auslieferung alle geltenden Vorschriften hinsichtlich elektromagnetischer Verträglichkeit (EMV) und Sicherheitsstandards.

Änderungen am Gerät können dazu führen, dass einer oder mehrere dieser EMV- und Sicherheitsstandards nicht mehr eingehalten werden. Änderungen an Ihrem Gerät beinhalten auch den Austausch von Teilen oder das Hinzufügen von Komponenten, Optionen oder Peripheriegeräten, die von Thermo Fisher Scientific nicht ausdrücklich für das Produkt autorisiert und freigegeben sind. Zur Sicherstellung der dauerhaften Einhaltung der EMV- und Sicherheitsstandards dürfen Ersatzteile und zusätzliche Komponenten, Zusatzmodule und Peripheriegeräte nur bei Thermo Fisher Scientific oder einer autorisierten Vertretung bestellt werden.

Das Gerät hat das Werk in einwandfreiem Zustand verlassen.

### Sehen Sie dazu auch

Informationen zur Konformität (> Seite 188)

# 3 Überblick über das Gerät

Dieses Kapitel stellt Ihnen die besonderen Merkmale des Geräts und die wichtigsten Komponenten vor.

# **3.1** Ausstattung des Detektors

Das Gerät ist im Wesentlichen wie folgt ausgestattet:

- Eine Xenon-Blitzlampe als Lichtquelle für den gesamten Anregungs-Wellenlängenbereich von 200 nm bis 880 nm
- Optional ein zweiter Photomultiplier (PMT) zur Erweiterung des Emissionswellenlängenbereichs in den nahen infraroten Bereich (bis zu 900 nm), ohne dass der Detektor an Empfindlichkeit im UV- /VIS-Bereich verliert
- Datenaufnahmeraten im Single-Channel-Modus bis zu 100 Hz (Detektoren vom Typ VC) bzw. 200 Hz (Detektoren vom Typ VF mit der Software Chromeleon 7)
- Wellenlängenschaltungen < 250 ms
- Höchste Empfindlichkeit (Signal-/Rauschverhältnis höher als 550 für das Raman-Spektrum von Wasser bei 350 nm Anregung, über die gesamte Lebensdauer der Lampe). Dadurch können auch kleinste Peaks erkannt werden.
- Messzellen mit aktiver Temperaturregelung, um eine verbesserte Reproduzierbarkeit bei schwankenden Umgebungstemperaturen zu gewährleisten
- Ein Langpassfilter zur Unterdrückung des bei Gitterspektrometern auftretenden Lichtanteils höherer Ordnungen und von Streulicht. Bei VF-Detektoren kann über einen Motor das Filterrad zwischen 5 Stellungen bewegt werden.
- Messungen von bis zu vier Daten-Kanälen gleichzeitig (nur VF-Detektoren) mit unabhängigen Parameter Einstellungen (PMT Auswahl, Wellenlängen, Sensitivity, Filterrad).

# 3.2 Funktionsprinzip

Fluoreszenz-Detektoren gehören zu den optischen Detektoren. In einem Fluoreszenz-Detektor wird die Probe mit Licht einer definierten Wellenlänge bestrahlt. Das Licht wird von der Probensubstanz absorbiert und die Substanz dadurch in einen angeregten Zustand versetzt (Anregung). Wenn die Probensubstanz ihren Grundzustand wieder annimmt, strahlt sie Licht mit einer größeren Wellenlänge ab (Emission). Der Fluoreszenz-Detektor erfasst das von fluoreszierenden Substanzen emittierte Licht. Dabei wird der Photomultipler (PMT) im 90°-Winkel zur Lichtquelle positioniert.

Im Gegensatz zu UV-/VIS-Detektoren wird also nicht die Differenz zwischen zwei Lichtintensitäten (Absorption) gemessen, sondern ein sehr schwaches Lichtsignal.



Abbildung 1: Vereinfachte Darstellung eines lichtabsorbierenden Moleküls



Abbildung 2: Vereinfachte Darstellung eines lichtemittierenden Moleküls (Fluoreszenz)

Beispiele für die Anwendung von Fluoreszenz im täglichen Leben sind Textmarker oder optische Aufheller in Waschmitteln ("Weißmacher"). Bei Textmarkern absorbieren Tagesleuchtfarben (Neonfarben) den blauen und nahen, nicht sichtbaren UV-Bereich des Tageslichtes und emittieren Licht größerer Wellenlänge (typisch sind blaugrün, gelb und rot).

Wie in der folgenden Abbildung gezeigt, wird der Lichtstrahl der Xenon-Blitzlampe (Nr. 1) über die Lampenoptik (Nr. 2) auf den Eingang des Anregungs-Monochromators (Nr. 3) gebündelt. Der Anregungs-Monochromator überträgt nur das Licht mit der ausgewählten Anregungs-Wellenlänge auf die Probe in der Messzelle (Nr. 4). Der Großteil des Lichts durchdringt die Probe und regt diese dabei zum Fluoreszieren an. Hinter der Messzelle misst ein Referenzsensor (Nr. 5) die Intensität des Anregungslichts. Das Referenzsignal dient zur Kompensation von Fluktuationen der Lampenintensität und verbessert so die Empfindlichkeit.

Nach dem Austritt aus der Messzelle wird das emittierte Licht über die Emissionsoptik (Nr. 6) auf den Emissions-Monochromator (Nr. 8) gebündelt.

Vor dem Emissions-Monochromator befindet sich ein Langpassfilter, welcher nur Licht oberhalb einer bestimmten Wellenlänge passieren lässt. In VF-Detektoren ist ein Filterrad (Nr. 7) installiert, das mit Hilfe eines Motors zwischen 5 Stellungen mit unterschiedlichen Grenzwellenlängen bewegt werden kann. Damit kann in vielen Anwendungen eine noch bessere Empfindlichkeit erreicht werden und zusätzlich erhöht sich die Flexibilität bei der Methodenentwicklung. VC-Detektoren besitzen einen Filter mit einer festen Grenzwellenlänge von 280 nm.

Der Emissions-Monochromator überträgt nur das Licht mit der ausgewählten Emissions-Wellenlänge auf einen Photomultiplier (PMT) (Nr. 9), wo die Messung der Lichtintensität erfolgt. Ein zweiter, (optionaler) rotsensitiver Photomultiplier (Nr. 10) misst Licht im nahen Infrarotbereich (bis 900 nm).



Abbildung 3: Aufbau	ı der Optik	(schematisch)
---------------------	-------------	---------------

Nr.	Komponente	Beschreibung
1	Xenon-Blitzlampe	Lichtquelle für die Wellenlängen vom UV-Bereich bis zum nahen Infrarot-Bereich
2	Lampenoptik	Bündelt den Lichtstrahl der Xenon-Blitzlampe auf den Eingang des Anregungs-Monochromators
3	Anregungs- Monochromator	Lässt nur Licht mit der ausgewählten Anregungs- Wellenlänge passieren
4	Messzelle	Der Eluent mit den Analyten fließt durch die Messzelle. Das Anregungslicht gelangt durch die Messzelle zum Referenzsensor, das Fluoreszenzlicht verlässt die Messzelle im 90°-Winkel dazu.
5	Referenzsensor	Misst das Anregungslicht durch die Messzelle und wird zur Kompensierung von Lampenfluktuationen genutzt
6	Emissionsoptik	Bündelt den emittierten Lichtstrahl von der Messzelle auf den Emissions-Monochromator
7	Filterrad	Trägt die optischen Filter zum Ausblenden von Licht unterhalb der Langpasswellenlänge des ausgewählten Filters
8	Emissions- Monochromator	Lässt nur Licht mit der ausgewählten Emissions- Wellenlänge passieren
9	Photomultiplier (PMT)	Wandelt Licht in einen messbaren Strom
10	Zweiter PMT	Misst Licht im nahen Infrarotbereich (bis 900 nm)

# 3.3 Innenansicht

Die für den Anwender zugänglichen Komponenten des Geräts befinden sich direkt hinter den Gerätetüren:



Abbildung 4: Innenansicht (hier mit eingebauter Messzelle)

Nr.	Beschreibung
1	Kühllufteinlass
2	Tastatur mit Statusanzeigen
3	Messzelle
4	Leckage-Wanne mit Leaksensor
5	Zwischenwand
	Die Aussparungen in der Zwischenwand werden mit Hilfe spezieller Einsätze zur Kapillarführung verwendet.
6	Typenschild mit Angabe von Modulnamen, Seriennummer, Teilenummer und Revisionsnummer (wenn vorhanden)
7	Leaksensor
8	Kapillarclips
# 3.4 Messzelle

Der Detektoraufbau ermöglicht es, leicht zur Messzelle im Innenraum des Detektors zu gelangen.



Abbildung 5: Messzelle (Beispiel)

Nr.	Beschreibung
1	Befestigungsschrauben der Messzelle - Zur Befestigung von der Messzelle an den Detektor.
2	Messzellen-Typenaufkleber
3	Auslass - Für den Anschluss der Waste-Verbindung.
4	Einlass - Für den Anschluss der Einlass-Kapillare.
5 5a, 5b, 5c	Optischer Block - Berühren Sie nie den optischen Block. Optische Ports

#### Messzellen-Typenaufkleber

An der Messzelle befinden sich ein oder mehrere Messzellen-Typenaufkleber, die Informationen über die Messzelle enthalten, wie zum Beispiel Messzellen-Typ, Artikelnummer und Seriennummer.

### Messzellen-Identifikationschip

Ein Identifikationschip (ID-Chip) an der Messzelle speichert Informationen, wie Messzellen-Typ und Seriennummer der Messzelle. Der ID-Chip speichert außerdem auch Daten während dem Betrieb, wie zum Beispiel die Zeit, die die Messzelle dem Lampenlicht ausgesetzt ist.

Wenn die Messzelle installiert ist und die Detektortüren geschlossen sind, liest das Gerät die Daten vom Chip und übermittelt die Messzellendaten an die Chromeleon-Software.

#### Temperaturreglung

Die Messzellen verfügen über eine Temperaturregelung. Messzelle und Wärmetauscher können auf eine benutzerdefinierte Temperatur erwärmt werden.

Der Wärmetauscher passt die Temperatur der mobilen Phase an die Temperatur der Messzelle an, ehe die mobile Phase in den optischen Flussweg innerhalb der Messzelle eintritt. Das hier angegebene Volumen von Wärmetauscher und Einlasskapillare beeinflusst die Retentionszeiten sowie Peakbreiten.

#### Messzellentypen

Alle Messzellen sind für schnelle Trennungen ohne Verlust der chromatographischen Auflösung optimiert. Beachten Sie auch die allgemeinen Hinweise zum Umgang mit Messzellen unter Hinweise zum Gebrauch von Messzellen () Seite 100).

Fragen zu den Messzellen oder zur Verfügbarkeit anderer Messzellen beantwortet Ihnen gern die Thermo Fisher Scientific-Vertriebsorganisation.

Bestellinformationen zu den Messzellen finden Sie unter Optionales Zubehör (> Seite 184).

Informationen zu den Messzellen-Spezifikationen wie auch zu den Materialien im Flussweg und den Temperaturbereich finden Sie unter Messzellen (> Seite 178).

# 3.5 Lampe

Als Lichtquelle dient eine Xenon-Blitzlampe.



Abbildung 6: Xenon-Blitzlampe

- Um die Lebensdauer der Lampe zu verlängern, wird diese erst automatisch eingeschaltet, wenn die Datenaufnahme beginnt, und nach Beenden der Datenaufnahme wieder ausgeschaltet.
- Die Lampe wird je nach gewähltem Lampen-Modus mit unterschiedlicher Blitzrate betrieben. Durch die Wahl eines anderen Lampenmodus in Phasen, in denen keine interessanten Peaks eluieren, kann die Lebensdauer der Lampe verlängert werden.
- Sie können die Betriebsdauer der Lampe überwachen. Diese Funktion soll helfen, festzustellen, wann eine Lampe ausgetauscht werden muss.
- Der Tausch der Lampe muss von einem Thermo Fisher Scientific-Servicetechniker durchgeführt werden.

Einzelheiten zu den verfügbaren Lampen-Modi und wie die Lebensdauer verlängert werden kann, finden Sie unter Lampen-Modus (> Seite 119).

Weitere Informationen zur Überwachung der Betriebsdauer der Lampe finden Sie unter Überwachen der Lampenbetriebsdauer () Seite 136).

# **3.6 Erkennen von Undichtigkeiten (Leakerkennung)**

Undichtigkeiten sind ein potentielles Sicherheitsrisiko.

Der Leaksensor im Gerät überwacht die fluidischen Verbindungen im Gerät auf Undichtigkeiten. Die Flüssigkeit sammelt sich in der Leckage-Wanne und wird zum Ablauf geleitet. Durch den Ablauf fließt die Flüssigkeit über das Drainagesystem des Vanquish-Systems in den Abfall.

Wenn der Leaksensor auf eine Undichtigkeit anspricht, leuchtet die Statusanzeige rot und ein akustisches Signal ertönt. Finden und beseitigen Sie die Ursache wie in diesem Handbuch beschrieben.

# 3.7 Betrieb

Das Gerät wird über einen Rechner gesteuert, auf dem das Chromatographie-Datensystem (CDS) Chromeleon installiert ist. Mit Hilfe der Software Chromeleon führen Sie die Gerätesteuerung und Datenaufnahme aus und verwalten Ihre Daten direkt in der Software.

Eine kurze Beschreibung der Gerätesteuerung und automatisierten Probenanalyse mit der Software Chromeleon finden Sie in der Vanquish-System-Betriebsanleitung. Einzelheiten zu Steuerbefehlen und Bedienung des Geräts finden Sie in der Chromeleon-Hilfe.

**TIPP** Das Gerät kann auch mit anderen Datensystemen, wie Xcalibur<sup>™</sup>, betrieben werden. Dabei ist die Installation weiterer Software zusätzlich zum Datensystem erforderlich. Weitere Fragen beantwortet Ihnen gern die Thermo Fisher Scientific-Vertriebsorganisation.

Im Gerät befindet sich eine Tastatur, mit der Sie direkt am Detektor das akustische Signal für den aktuellen Alarm abschalten sowie den Detektor initialisieren können.

# 4 Auspacken

In diesem Kapitel finden Sie Informationen rund um das Auspacken des Geräts sowie zum Lieferumfang.

# 4.1 Auspacken

#### Beschädigte Verpackung, Mängel bei Ankunft des Geräts

Überprüfen Sie die Transportverpackung auf Anzeichen äußerer Beschädigung und überprüfen Sie das Gerät nach dem Auspacken auf Anzeichen mechanischer Beschädigungen, die auf dem Versandweg aufgetreten sein könnten.

Besteht der Verdacht, dass das Gerät auf dem Versandweg in irgendeiner Weise beschädigt wurde, melden Sie etwaige Schäden sofort sowohl dem Transportunternehmen als auch Thermo Fisher Scientific. Nur bei sofortiger Reklamation kommt die Transportversicherung für die aufgetretenen Schäden auf.

#### Auspacken des Geräts



#### VORSICHT—Schweres und unhandliches Gerät

Das Gerät ist zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden am Gerät zu vermeiden, beachten Sie folgende Hinweise:

- Um das Gerät zu greifen und zu transportieren, sind mindestens zwei Personen erforderlich, zum Beispiel, um das Gerät anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn das Gerät in den Systemturm gehoben oder aus dem Systemturm entfernt wird.
- Verwenden Sie f
  ür den Transport des Ger
  äts die Tragegriffe, die mitgeliefert wurden. Transportieren Sie oder heben Sie das Ger
  ät niemals an den Ger
  ätet
  üren. Dies f
  ührt zu Besch
  ädigungen an den T
  üren oder am Ger
  ät.

#### Erforderliche Werkzeuge

Schraubendreher, Torx<sup>™</sup> T20

Gehen Sie wie folgt vor

- 1. Stellen Sie den Versandkarton auf den Boden und öffnen Sie ihn.
- 2. Entnehmen Sie das Zubehör.
- Nehmen Sie das Gerät aus dem Versandkarton heraus: Greifen Sie das Gerät dazu an den Tragegriffen. Heben Sie das Gerät langsam und vorsichtig aus dem Versandkarton heraus.



Abbildung 7: Tragegriffe am Gerät

Nr.	Komponente
1	Tragegriffe
2	Befestigungsschraube (eine Schraube an jedem Tragegriff)

- 4. Stellen Sie das Gerät auf eine stabile Oberfläche.
- Wenn zutreffend: Entfernen Sie sonstiges Verpackungsmaterial. Entfernen Sie etwaige Schutzfolien erst dann von den Oberflächen des Geräts, wenn dieses korrekt im Systemturm positioniert ist.
- Greifen Sie das Gerät an den Tragegriffen und transportieren Sie es zum Aufstellungsort, falls es sich noch nicht dort befindet, und positionieren Sie das Gerät im Systemturm (siehe Systemaufbau (▶ Seite 54)).
- 7. Lösen Sie die Befestigungsschraube an jedem der beiden Tragegriffe, bis Sie den Tragegriff in der Schiene bewegen können. Entfernen Sie die Schrauben nicht vollständig von den Tragegriffen.
- 8. Ziehen Sie die Tragegriffe zur Rückseite des Geräts hin aus den Schienen heraus.



Abbildung 8: Herausziehen des Tragegriffs aus der linken Schiene

**TIPP** Bewahren Sie den Versandkarton, die Tragegriffe mit den Befestigungsschrauben und die Verpackungsmaterialien auf. Sie benötigen diese Dinge, wenn Sie das Gerät an einen anderen Ort transportieren oder verschicken möchten.

9. Einige Oberflächen wie die Türen des Geräts sind für den Transport durch Folien geschützt. Entfernen Sie die Schutzfolien soweit vorhanden von allen Oberflächen.

# 4.2 Lieferumfang

Im Lieferumfang enthalten sind:

- Detektor
- Zubehörkit
- Betriebsanleitung
- Netzkabel

Informationen zum Inhalt des Zubehörkits sowie zur Nachbestellung finden Sie unter Zubehör, Ersatzteile und Verbrauchsmaterialien (▶ Seite 181).

# **5** Installation

In diesem Kapitel finden Sie Informationen zu den Anforderungen an den Aufstellungsort sowie zum Aufbau, zur Installation und zur Konfiguration des Geräts im Vanquish-System und in der Chromatographie-Software.

# 5.1 Sicherheitshinweise für die Installation

Beachten Sie folgende Sicherheitshinweise:



Beachten Sie alle Warnhinweise und Vorsichtsmaßnahmen in Sicherheitsmaßnahmen (▶ Seite 21).



#### VORSICHT—Schweres und unhandliches Gerät

Das Gerät ist zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden am Gerät zu vermeiden, beachten Sie folgende Hinweise:

- Um das Gerät zu greifen und zu transportieren, sind mindestens zwei Personen erforderlich, zum Beispiel, um das Gerät anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn das Gerät in den Systemturm gehoben oder aus dem Systemturm entfernt wird.
- Verwenden Sie f
  ür den Transport des Ger
  äts die Tragegriffe, die mitgeliefert wurden. Transportieren Sie oder heben Sie das Ger
  ät niemals an den Ger
  ätet
  üren. Dies f
  ührt zu Besch
  ädigungen an den T
  üren oder am Ger
  ät.



#### VORSICHT—Stromschlag oder Schäden am Gerät

Auch im ausgeschalteten Zustand fließt im Gerät Strom, solange das Netzkabel eingesteckt ist. Reparaturen am Gerät, während das Gerät am Stromnetz angeschlossen ist, können zu Personenschäden führen.

- Ziehen Sie immer das Netzkabel ab, bevor Sie Reparaturen im Geräteinneren durchführen.
- Sollten Sie Gehäuseabdeckungen oder Seitenwände entfernen müssen, schließen Sie das Netzkabel keinesfalls an das Gerät an, solange die Abdeckungen und Seitenwände noch nicht montiert sind.

# 5.2 Installieren des Gerätes

Ein Servicetechniker von Thermo Fisher Scientific installiert das Vanquish-System einschließlich aller mitgelieferten Module, Optionen und Teile, und nimmt das System in Betrieb. Der Servicetechniker prüft, dass das Vanquish-System korrekt installiert wurde und alle Module sowie das gesamte System unter Einhaltung der Spezifikationen funktionieren. Der Servicetechniker führt zudem die Grundfunktionen und wichtigsten Merkmale vor.

Wenn der Aufbau des Geräts durch eine andere Person als einen Thermo Fisher Scientific-Servicetechniker erfolgt, folgen Sie den Schritten in diesem Kapitel.

#### ACHTUNG

Das Gerät ist Teil des Vanquish-Systems. Befolgen Sie daher die Reihenfolge für die Installation der Systemmodule, die in der *Betriebsanleitung zum Vanquish-System* beschrieben ist.

- Beachten Sie die Sicherheitshinweise und Anforderungen an den Aufstellungsort. Siehe Sicherheitshinweise für die Installation (▶ Seite 48) und Anforderungen an den Aufstellungsort (▶ Seite 51).
- 2. Bauen Sie das Gerät auf. Siehe Aufbauen der Hardware (> Seite 54).
- 3. Stellen Sie die Flussverbindungen her. Siehe Anschließen der Flussverbindungen (▶ Seite 64).
- Schalten Sie das Gerät ein. Siehe Einschalten des Gerätes (▶ Seite 88).

#### TIPP

Bevor Sie ein Modul des Vanquish-Systems zum ersten Mal einschalten vergewissern Sie sich, dass die Chromatographie-Software auf dem Datensystemrechner installiert ist. Die erforderlichen USB-Treiber werden automatisch geladen und das Windows™-Betriebssystem kann das Gerät erkennen, wenn dieses eingeschaltet ist.

- 5. Setzen Sie das Gerät in der Software auf. Siehe Einrichten des Gerätes in der Software (▶ Seite 89).
- 6. Führen Sie eine Wellenlängenkalibrierung und eine Wellenlängenvalidierung durch.

7. Empfohlen:

Führen Sie eine Qualifizierung über Instrument Installation Qualification durch.

In der Software Chromeleon führt Sie ein Assistent durch den Qualifizierungsvorgang. In der **Chromeleon 7 Console**: Klicken Sie auf **Tools > Instrument Qualification > Installation Qualification**.

Folgen Sie den Anweisungen in der Bedienungsanleitung zur Instruments Installation Qualification. Die Anleitung enthält alle Informationen zum benötigten Material sowie detaillierte Anweisungen.

#### ACHTUNG

Wird das Gerät mit einem anderen Datensystem betrieben, lesen Sie in der Dokumentation zu der verwendeten Software nach und/oder führen Sie die Qualifizierung manuell durch. Die *Bedienungsanleitung zur Instruments Installation Qualification* enthält Informationen zu den Parametern, die angepasst werden müssen, sowie zu den erforderlichen Einstellungen.

 Empfohlen: Führen Sie eine Qualifizierung über Operational Qualification durch.
 Das Qualifizierungskit enthält alle erforderlichen Materialien für die Qualifizierung sowie detaillierte Anweisungen.

#### Transportieren des Geräts nach der Installation

Falls Sie das Gerät transportieren müssen, nachdem es aufgestellt und im Vanquish-System installiert wurde, bereiten Sie es für den Transport vor und transportieren Sie es an den neuen Aufstellungsort. Folgen Sie den Anweisungen in Transportieren und Versenden des Gerätes () Seite 152).

#### Sehen Sie dazu auch

- Durchführung einer Wellenlängen-Kalibrierung (> Seite 137)
- Durchführung einer Wellenlängenvalidierung (> Seite 139)

# 5.3 Anforderungen an den Aufstellungsort

Die Umgebungsbedingungen sind wichtig, um den optimalen Betrieb des Gerätes zu ermöglichen. Dieser Abschnitt behandelt wichtige Anforderungen an den Aufstellungsort. Beachten Sie Folgendes:

- Betreiben Sie das Gerät nur bei angemessenen Laborbedingungen.
- Das Gerät ist zur Verwendung im Vanquish-System gedacht. Beachten Sie die Anforderungen an den Aufstellungsort für das Vanquish-System in der *Vanquish-System-Betriebsanleitung*.
- Die Spezifikationen finden Sie unter Spezifikationen (▶ Seite 175) und im Abschnitt *Spezifikationen* in den *Betriebsanleitungen* für die anderen Module im Vanquish-System.
- Zu den allgemeinen Restrisiken, siehe Allgemeine Restrisiken (▶ Seite 24).

### 5.3.1 Hinweise zur Stromversorgung

Das Netzteil des Geräts verfügt über einen weiten Eingangsspannungsbereich und kann sich automatisch an jede Netzspannung innerhalb des spezifizierten Bereichs für das Gerät anpassen.



#### VORSICHT—Stromschlag oder Schäden am Gerät

Wird das Gerät an höhere oder niedrigere als die angegebenen Spannungen angeschlossen, kann dies zu Personenschäden oder Schäden am Gerät führen.

Schließen Sie das Gerät nur an die angegebene Netzspannung an.

### 5.3.2 Netzkabel

Die Netzkabel sind den länderweise unterschiedlichen Wandsteckdosen angepasst. Die Buchse, die an den Netzstecker des Geräts angeschlossen wird, ist bei allen Netzkabeln gleich. Der Stecker des Netzkabels, der an die Wandsteckdose angeschlossen wird, ist unterschiedlich.



#### WARNUNG—Stromschlag oder Schäden am Gerät

- Verwenden Sie nur die von Thermo Fisher Scientific für das Gerät bereitgestellten Netzkabel.
- Verwenden Sie ausschließlich ein Netzkabel, das für das Land bereitgestellt wurde, in dem Sie das Gerät betreiben.
- Verwenden Sie keine Verlängerungskabel.
- Schließen Sie das Netzkabel niemals an eine Steckdose an, an die auch andere Geräte angeschlossen sind (zum Beispiel Mehrfachsteckdosen).
- Betreiben Sie Ihr Gerät nur an einer Spannungsquelle mit Schutzerdung.
- Im Notfall muss das Netzkabel des Gerätes einfach zugänglich sein, damit Sie das Gerät jederzeit vom Stromnetz trennen können.



#### WARNUNG—Stromschlag oder Schäden an einem Produkt

Unzweckmäßiger Gebrauch von Netzkabeln kann zur Gefährdung Ihrer Person oder Schäden am Gerät führen. Verwenden Sie die Netzkabel von Thermo Fisher Scientific ausschließlich für den Zweck, für den sie bestimmt sind. Verwenden Sie die Netzkabel nicht für andere Zwecke, zum Beispiel das Anschließen von anderen Geräten.

#### 5.3.3 Kondensation

# ACHTUNG—Kondensation im Geräteinneren kann die Elektronik und die Optik beschädigen.

- Vermeiden oder minimieren Sie bei Betrieb, Versand oder Lagerung Bedingungen, die zu einer Kondensatbildung im Gerät führen können. Vermeiden Sie, zum Beispiel, signifikante und schnelle Veränderungen der Umgebungsbedingungen.
- Besteht der Verdacht, dass sich Kondenswasser gebildet hat, lassen Sie das Gerät akklimatisieren. Dies kann einige Stunden dauern.
   Warten Sie, bis sich das Kondenswasser vollständig verflüchtigt hat, bevor Sie das Gerät an das Stromnetz anschließen.

# 5.4 Öffnen des Innenraums

Um zu den Komponenten im Innenraum des Gerätes zu gelangen, öffnen Sie die Gerätetüren. Direkt hinter den Gerätetüren sind die für den Anwender zugänglichen Komponenten und Flussverbindungen des Gerätes leicht erreichbar.



Abbildung 9: Öffnen der Gerätetüren

# 5.5 Aufbauen der Hardware

In diesem Abschnitt finden Sie Informationen, wie die Hardware aufgebaut werden muss, und welche Anschlüsse und Kabel für das Gerät notwendig sind.

### 5.5.1 Systemaufbau

Das Gerät ist Teil des Vanquish-Systems. Die Module des Systems werden typischerweise in einem Systemturm aufgebaut, wobei die genaue Anordnung von der Systemkonfiguration abhängt.

Die unten aufgeführten Abbildungen zeigen Konfigurationen mit einem einzelnen Fluoreszenz-Detektor sowie mit einem Fluoreszenz-Detektor als zweiten Detektor oberhalb eines UV-/VIS-Detektors.

#### TIPP

Da die Messzelle nur einen geringen Gegendruck aushält, sollte der Fluoreszenz-Detektor wann immer möglich das letzte Gerät im Flussweg sein.

Informationen zum Aufbau des Systems können Sie der Vanquish-System-Betriebsanleitung entnehmen.



System mit einem einzelnen Detektor



Nr.	Beschreibung
1	Solvent Rack
2	Fluoreszenz-Detektor
3	Autosampler
4	Pumpe
5	Systemsockel
6	Säulenthermostat



System mit einem Fluoreszenz-Detektor als zweiten Detektor



Nr.	Beschreibung
1	Solvent Rack
2	Fluoreszenz-Detektor
3	UV-/VIS-Detektor
4	Autosampler
5	Pumpe
6	Systemsockel
7	Säulenthermostat

## 5.5.2 Verbinden des Gerätes

#### Anschlüsse am Gerät

Folgende Anschlüsse befinden sich am Gerät:





Nr.	Beschreibung
1	Typenschild mit Angabe von Seriennummer, Teilenummer, Modulnamen, Revisionsnummer (wenn vorhanden), Netzspannung und Nennleistung sowie Herstelleradresse
2	Hauptnetzschalter (Ein/Aus)
3	Sicherungshalter
4	Netzbuchse
5	System Interlink-Port Für das Ein- und Ausschalten des Detektors über den Vanquish- Systemsockel sowie die Kommunikation bzw. Synchronisation zwischen dem Detektor und anderen Modulen im Vanquish-System. Zum Beispiel ermöglicht die Verbindung zwischen Autosampler und Detektor eine automatische Synchronisation der Probeninjektion mit dem Beginn der Datenaufnahme im Detektor. Die Synchronisation verbessert dadurch die Reproduzierbarkeit der Retentionszeit.

Nr.	Beschreibung
6	Digital I/O-Ports (Dig I/O)Für den Austausch digitaler Signale mit externen Geräten
	Jeder digitale E/A-Port verfügt über einen digitalen Eingang und einen Relaisausgang. Informationen zum Anschluss und zur Steckerbelegung finden Sie in Digital I/O (▶ Seite 191).
7	USB-Hub (Stecker Typ "A") Für den Anschluss anderer Module des Vanquish-Systems
8	USB-Port (Universal Serial Bus) (Stecker Typ "B") Für den Anschluss an andere Module im Vanquish-System oder an den Rechner, auf dem das Datenmanagement-System, zum Beispiel die Software Chromeleon, installiert ist

**TIPP** Thermo Fisher Scientific empfiehlt, die USB-Ports nur wie oben beschrieben zu verwenden. Werden die USB-Ports für andere Zwecke verwendet, kann Thermo Fisher Scientific den ordnungsgemäßen Betrieb nicht gewährleisten.

#### Gehen Sie wie folgt vor

#### ACHTUNG

- Verwenden Sie keine beschädigten Kommunikationskabel. Tauschen Sie das Kabel aus, wenn Sie den Verdacht haben, das Kabel sei beschädigt.
- Verwenden Sie zum Anschluss des Geräts ausschließlich die von Thermo Fisher Scientific zur Verfügung gestellten Kabel, um eine einwandfreie Funktion der Verbindung zu gewährleisten.
- 1. Platzieren Sie das Gerät so im System, wie dies entsprechend der Systemkonfiguration vorgegeben ist. Einzelheiten hierzu finden Sie in der *Vanquish-System-Betriebsanleitung*.
- 2. Schließen Sie die erforderlichen Verbindungskabel an das Gerät an.
- Schließen Sie das Netzkabel an (siehe Anschließen des Netzkabels (▶ Seite 61)).

#### Sehen Sie dazu auch

Anschließen der Verbindungskabel (> Seite 59)

### 5.5.2.1 Anschließen der Verbindungskabel

Der Anschluss der Verbindungskabel hängt davon ab, ob der Detektor der einzige Detektor im Vanquish-System ist oder der zweite.

#### Der Detektor ist der einzige Detektor im System

Schließen Sie die erforderlichen Verbindungskabel an den Detektor an. Informationen dazu, wie Sie den Detektor mit den anderen Modulen im Vanquish-System oder mit dem Chromatographie-Datensystemrechner verbinden, finden Sie in der *Vanquish-System-Betriebsanleitung*.

#### Der Detektor ist der zweite Detektor im System

Wenn der Fluoreszenz-Detektor der zweite Detektor im Vanquish-System ist (zum Beispiel nach dem Diodenarray-Detektor), stellen Sie die USB- und System-Interlink-Verbindungen entsprechend der Abbildung her.

- 1. Folgen Sie den Anweisungen zum Verbinden der anderen Module einschließlich des Diodenarray-Detektors in der *Vanquish-System-Betriebsanleitung*.
- 2. Verbinden Sie ein USB-Kabel von einem freien USB-Port am Diodenarray-Detektor mit dem Fluoreszenz-Detektor.
- Verbinden Sie ein System-Interlink-Kabel vom freien System Interlink-Port am Diodenarray-Detektor mit dem Fluoreszenz-Detektor.



4. Verbinden Sie ein System-Interlink-Kabel vom freien **System** Interlink-Port am Diodenarray-Detektor mit dem Säulenthermostat.

Abbildung 13: Kabelverbindungen im Vanquish-System mit Fluoreszenz-Detektor und Diodenarray-Detektor (Beispiel)

Nr.	Beschreibung
1	Systemsockel
2	Pumpe
3	Autosampler
4	Diodenarray-Detektor
5	Fluoreszenz-Detektor
6	Säulenthermostat
7	Verbindung zum Rechner

## 5.5.2.2 Anschließen des Netzkabels

### ACHTUNG

Kondensation im Gerät kann die Elektronik beschädigen.

- Vergewissern Sie sich, dass sich das Kondenswasser vollständig verflüchtigt hat, bevor Sie die Geräte an das Stromnetz anschließen.
- Besteht der Verdacht, dass sich Kondenswasser gebildet hat, lassen Sie das Gerät langsam akklimatisieren. Warten Sie, bis sich das Kondenswasser vollständig verflüchtigt hat, ehe Sie fortfahren.
- 1. Vergewissern Sie sich, dass alle Geräte über den Netzschalter ausgeschaltet sind.
- 2. Verbinden Sie das Netzkabel mit der Netzbuchse am Gerät.
- 3. Verbinden Sie das andere Ende des Netzkabels mit einer geeigneten Stromquelle.

### 5.5.3 Einbauen der Messzelle

Dieser Abschnitt beschreibt die Installation der Messzelle im Zuge der Erstinstallation des Detektors.

Anweisungen zum Ausbau einer Messzelle oder der Installation einer Messzelle nach der Aufbewahrung siehe Messzelle (> Seite 141).

#### ACHTUNG

Die Messzellen reagieren äußerst empfindlich auf Verschmutzung und Staub. Beachten Sie die folgenden Hinweise beim Einbau der Messzelle in den Detektor:

- Berühren Sie nie den optischen Block der Messzelle oder die empfindliche Elektronik auf der Rückseite.
- Die optischen Ports an der Messzelle sind empfindlich gegen Verunreinigungen und Kratzern. Berühren Sie niemals die optischen Ports der Messzelle und tauchen Sie diese niemals unter.
- Um Schäden an den optischen Ports der Messzelle zu vermeiden, gehen Sie vorsichtig vor, wenn Sie die Messzelle in den Messzellenschacht des Geräts einschieben.
- Verwenden Sie die vorgesehene Verpackung für die Messzelle, wenn Sie sie aufbewahren oder transportieren.
- Auf der Rückseite der Messzelle befinden sich die empfindlichen Kontaktpunkte für den Identifikationschip. Berühren Sie niemals die Kontaktpunkte. Vermeiden Sie Schäden an der Elektronik des ID-Chips.

#### Erforderliche Teile

Messzelle

#### Vorbereitungen

1. Entfernen Sie die Abdeckung vom Messzellenschacht. Lösen Sie dazu die beiden Schrauben per Hand. Die Schrauben sind unverlierbar mit der Abdeckung verbunden und müssen nicht entfernt werden.

**TIPP** Bewahren Sie die Abdeckung auf, um den Messzellenschacht zu verschließen, wenn keine Messzelle im Gerät eingebaut ist, insbesondere zum Transport oder Versand des Detektors.

2. Packen Sie die Messzelle aus.





Nr.	Beschreibung
1	Befestigungsschrauben der Messzelle - Zur Befestigung von der Messzelle an den Detektor.
2	Messzellen-Typenaufkleber
3	Auslass - Für den Anschluss der Waste-Verbindung.
4	Einlass - Für den Anschluss der Einlass-Kapillare.
5	Optischer Block - Berühren Sie nie den optischen Block.
5a, 5b, 5c	Optische Ports

#### Gehen Sie wie folgt vor

1. Setzen Sie die Messzelle gerade in den Messzellenschacht ein.



Abbildung 15: Einsetzen der Messzelle

2. Ziehen Sie die Befestigungsschrauben der Messzelle handfest an.

# 5.6 Anschließen der Flussverbindungen

### 5.6.1 Allgemeine Informationen und Hinweise

Beachten Sie folgende Regeln und Empfehlungen, wenn Sie Flussverbindungen anschließen:



Flussverbindungen können mit gefährlichen Substanzen gefüllt sein. Beachten Sie die Warnhinweise und Vorsichtsmaßnahmen in Sicherheitsmaßnahmen (> Seite 21).

#### ACHTUNG

Schmutzpartikel von anderen Modulen und Komponenten im System können sich in der Messzelle ablagern und sie verstopfen.

- Bevor Sie die Flussverbindungen zur Messzelle herstellen, f
  ühren Sie eine Kapillare vom System-Auslass in den Abfall und sp
  ülen Sie das System ohne den Detektor.
- Wenn Sie Geräte oder Komponenten im System installieren, spülen Sie diese immer in den Abfall, bevor Sie diese in den Flussweg des Systems aufnehmen. Folgen Sie den Anweisungen zum Spülen der Vanquish-Module in der Vanquish-System-Betriebsanleitung.

#### ACHTUNG

Messzellen reagieren äußerst empfindlich auf Verunreinigungen, Verstopfung und hohe Rückdrücke. Schon eine kurzzeitige Überschreitung des Maximaldrucks kann die Messzelle dauerhaft beschädigen. Beachten Sie die folgenden Hinweise, wenn Sie die Messzelle im System-Flussweg anschließen:

- Wenn Sie eine Komponente im Flussweg nach der Messzelle anschließen, beachten Sie den für die Messzelle spezifizierten Rückdruck.
- Verwenden Sie nur saubere Viper-Kapillaren, die mit der Messzelle mitgeliefert wurden und zuvor durch eine Verschlusskappe ordentlich verschlossen waren.
- Verwenden Sie ausschließlich die für die Messzelle bereitgestellte Waste-Leitung.
- Vermeiden Sie ein Verstopfen der Messzelle oder der Waste-Verbindung.

#### Gehen Sie wie folgt vor

Folgen Sie diesen Schritten, um die Flussverbindungen herzustellen und die Installation des Geräts abzuschließen:

- 1. Stellen Sie die Flussverbindungen her (siehe Flussverbindungen an der Messzelle (▶ Seite 71)).
- 2. Verbinden Sie das Gerät mit dem Drainage-System (Informationen hierzu finden Sie in der *Vanquish-System-Betriebsanleitung*).

Eine detaillierte Anleitung für die Installation und Hinweise zum Umgang finden Sie unter Anschließen von Fittingen, Kapillaren und Schläuchen (▶ Seite 68).

### 5.6.2 Kapillar- und Schlauchführung durch das System

Flussverbindungen zwischen den Modulen im Vanquish-System werden entweder durch den Schlauchkanal in den Geräten oder die Führungslöcher oder Kapillarclips der Geräte geführt.

#### Schlauchkanal, Schlauchführung, Schlauchklammer

Um Schläuche und Verbindungen (Schläuche für Lösungsmittel und Waschflüssigkeiten, Detektor-Waste-Verbindung) vom obersten Modul zum untersten Modul durch den Vanquish-Systemturm zu führen, sind die stapelfähigen Module innen rechts mit einem Schlauchkanal ausgestattet.

Der Schlauchkanal hat vier Schlauchführungen. Jede Führung kann bis zu drei Schläuche oder Verbindungen aufnehmen.

Drücken Sie den Schlauch (oder die Verbindung) in jedem Modul in die entsprechende Führung.



Abbildung 16: Schlauchkanal mit Schlauchführungen (links: Ansicht von innen, rechts: von oben)

Nr.	Zur Verwendung für
1	Ansaugschläuche für Lösungsmittel (bis zu drei Schläuche)
2	Ansaugschläuche für Lösungsmittel (bis zu drei Schläuche)
3	Schläuche für die Waschflüssigkeiten (Hinterspülung, Waschen der Autosampler-Nadel)
4	Detektor-Waste-Verbindung

Es stehen Schlauchklammern zur Verfügung, um die Schläuche in ihrer Position zu halten. Schieben Sie die Seite mit der Klammer auf die Drainage-Leitung.



Abbildung 17: Schlauchklammer (links); Schlauchklammer installiert (rechts)

Duale Systemaufbauten

Bei Dual-Systemaufbauten kann die Anzahl der Schläuche die Fassungskapazität der Schlauchführungen übersteigen. In diesem Fall wird empfohlen, die Lösungsmittelschläuche in den Schlauchführungen zu platzieren und andere Schläuche frei im Schlauchkanal entlang zu führen.

#### Führungslöcher und Kapillarclips

Die Systemmodule haben an bestimmten Positionen Führungslöcher und Kapillarclips. Führen Sie Flussverbindungen von einem Modul im Vanquish-System zum nächsten Modul durch das entsprechende Führungsloch oder den Kapillarclip, wenn Sie in dieser Anleitung dazu aufgefordert werden.

### 5.6.3 Installieren der Einsätze für die Zwischenwand

Im Zubehörkit des Detektors stehen zwei Einsätze für die Zwischenwand zur Verfügung.



Abbildung 18: Verfügbare Einsätze für die Zwischenwand

Nr.	Beschreibung
1	Einsatz mit Schlitz für Kapillaren mit einem kleinen Außendurchmesser, wie unisolierte Kapillaren.
2	Dreheinsatz für Kapillaren, die nicht in den Einsatz mit Schlitz passen, wie isolierte Kapillaren.

Installieren des Einsatzes mit Schlitz

- 1. Drücken Sie den Einsatz mit Schlitz in die erforderliche Aussparung der Zwischenwand (wenn nicht bereits vorinstalliert).
- 2. Stecken Sie zur sicheren Befestigung die Nase in die vorgesehene kleine Öffnung in der Zwischenwand.



Abbildung 19: Befestigung des Einsatzes mit Schlitz



Abbildung 20: Kapillare im Einsatz mit Schlitz

Installieren des Dreheinsatzes

- 1. Drücken Sie den den Dreheinsatz in die erforderliche Aussparung der Zwischenwand (wenn nicht bereits vorinstalliert).
- 2. Drehen Sie den Dreheinsatz nach vorn, um den Einsatz zu öffnen und anschließend die Kapillare durch den Einsatz zu führen.
- 3. Drehen Sie den Dreheinsatz Richtung Zwischenwand und schließen Sie so die Öffnung des Dreheinsatzes, damit die Kapillare nicht herausrutschen kann.



Abbildung 21: Verwendung des Dreheinsatzes

#### 5.6.4 Anschließen von Fittingen, Kapillaren und Schläuchen

Die Einlass- und Auslassports der Messzellen unterstützen die folgenden Kapillarfittinge:

- Viper™
- nanoViper™
- Standard 1/16" HPLC-Fittinge mit Schneidringen (PEEK; SST) oder fingerfeste Fittinge (PEEK)

In diesem Abschnitt finden Sie Informationen zum Anschluss von und Umgang mit Kapillaren, Fittingen und Schläuchen.

### 5.6.4.1 Allgemeine Hinweise

Folgen Sie beim Anschließen von Kapillaren und Schläuchen diesen allgemeinen Empfehlungen:

- Verwenden Sie nur die Kapillaren und Schläuche (zum Beispiel Lösungsmittelschläuche oder Ablaufschläuche), die mit dem Produkt mitgeliefert werden, oder die von Thermo Fisher Scientific als Ersatzteile oder optionales Zubehör empfohlen werden.
- An den Anschlüssen dürfen keine Verunreinigungen haften.
   Eindringende Schmutzpartikel können zu Schäden am System oder falschen Testergebnissen führen.
- Verwenden Sie keine übermäßig beanspruchten, eingekerbten, geknickten oder anderweitig beschädigten Kapillaren oder Schläuche.
- Installieren Sie Kapillarverbindungen nur an den dafür vorgesehenen Positionen.

### 5.6.4.2 Anschließen von Viper-Kapillaren

Dieser Abschnitt beschreibt den Anschluss von Viper™-Kapillaren. Alle Viper-Flussverbindungen im Vanquish-System sind so konzipiert, dass sie mit den Fingern ohne Werkzeug angezogen werden können.

Gehen Sie wie folgt vor, um Viper-Kapillaren mit Rändelschraube anzuschließen:

#### ACHTUNG

- Lösen und ziehen Sie die Viper-Kapillaren nur mit den Fingern fest. Verwenden Sie kein Werkzeug außer der Rändelschraube, die mit der Kapillare mitgeliefert wird.
- Um Schäden an den Kapillaren und Anschlüssen zu vermeiden, lösen und ziehen Sie die Viper-Kapillaren nur fest, wenn der Systemdruck gleich Null ist.



#### Abbildung 22: Viper-Fitting mit Rändelschraube

Nr.	Beschreibung
1	Rändelschraube
2	Kapillare
3	Schlitz

- 1. Führen Sie die Viper-Kapillare in den Anschlussport ein.
- 2. Ziehen Sie die Verbindung mit Hilfe der Rändelschraube fest.

**TIPP** Achten Sie auf den Schlitz in der Rändelschraube. Durch diesen Schlitz können Sie bei engen Verbindungen die Rändelschraube von benachbarten Verbindungen leicht entfernen und später wieder aufsetzen.

 Prüfen Sie die Verbindung auf Undichtigkeit. Wenn eine Undichtigkeit vorhanden ist, folgen Sie den Schritten weiter unten.

#### Beheben von Undichtigkeiten bei Viper-Fittingen mit Rändelschraube

- 1. Ziehen Sie die Verbindung etwas fester an.
- 2. Bleibt die Undichtigkeit bestehen, entfernen Sie die Kapillare.
- 3. Reinigen Sie die Kapillarenden vorsichtig mit einem fusselfreien Tuch, das mit Isopropanol angefeuchtet ist.
- 4. Schließen Sie die Kapillare wieder an.
- 5. Verwenden Sie eine neue Viper-Kapillare, wenn die Undichtigkeit weiter bestehen bleibt.

### 5.6.5 Flussverbindungen an der Messzelle

Verbinden Sie die Einlasskapillare und die Waste-Verbindung mit der Messzelle, wenn die Messzelle im Gerät eingebaut ist.

#### ACHTUNG

Rückdrücke, die über dem spezifizierten, maximalen Grenzwert der Messzelle liegen, können die Messzelle zerstören. Beachten Sie folgende Hinweise:

- Verwenden Sie ausschließlich die Waste-Leitung für Ihren Detektor.
- Stellen Sie die Waste-Verbindung zur Messzelle nur wie in der Anleitung beschrieben her.
- Leiten Sie keine Abfall-Flüssigkeiten von der Messzelle in den offenen Ablauf der Systemdrainage im Vanquish-System.
- Setzen Sie die Messzelle niemals einem übermäßigen Rückdruck aus.
- Vermeiden Sie ein Verstopfen der Messzelle oder der Waste-Verbindung.
- Wenn Sie eine Komponente im Flussweg nach der Messzelle anschließen, beachten Sie den für die Messzelle spezifizierten Rückdruck.

Beachten Sie zusätzlich dazu die Hinweise zum ordnungsgemäßen Anschluss der Messzelle in Allgemeine Informationen und Hinweise (> Seite 64).

#### Erforderliche Teile

- Einlasskapillare, abhängig von Ihrem Systemaufbau:
  - Der Detektor ist der einzige Detektor im Vanquish-System
     Verwenden Sie die Einlasskapillare vom System-Zubehörkit.
  - Der Detektor ist der zweite Detektor im Vanquish-System
     Verwenden Sie die mit dem Detektor mitgelieferte Kapillare mit größerem Innendurchmesser.
- Detektor-Waste-Leitung
   Folgen Sie den Anweisungen zum Herstellen der Waste-Verbindung unter Anschließen der Detektor-Waste-Leitung (> Seite 76).

#### Erforderliche Werkzeuge

Für die Detektor-Waste-Leitung: Kapillarschneider (optional)

Vorbereitungen

- 1. Führen Sie die Kapillaren von den Modulen im System-Flussweg vor dem Gerät in den Abfall und spülen Sie das System ohne das Gerät, bevor Sie die Flussverbindungen zur Messzelle herstellen. Informieren Sie sich in der *Vanquish-System-Betriebsanleitung*.
- 2. Entfernen Sie die Verschlussstopfen vom Einlass und Auslass der Messzelle.

**TIPP** Bewahren Sie die Verschlussstopfen der Messzelle auf, beispielsweise in der Messzellen-Verpackung, um sie für die Aufbewahrung oder den Transport der Messzelle zur Hand zu haben.

 Installieren Sie die Einsätze für die Zwischenwand (wenn sie noch nicht installiert wurden). Siehe Installieren der Einsätze für die Zwischenwand (▶ Seite 67).

#### Gehen Sie wie folgt vor

- 1. Schließen Sie die Einlasskapillare an die Messzelle an.
- 2. Schließen Sie die Detektor-Waste-Leitung an den Detektor an.

#### Sehen Sie dazu auch

- Anschließen der Einlasskapillare (> Seite 72)
- Anschließen der Detektor-Waste-Leitung (> Seite 76)

#### 5.6.5.1 Anschließen der Einlasskapillare

Abhängig von den Modulen in Ihrem Vanquish-System können Sie die Einlasskapillare direkt vom Säulenthermostat oder von einem vorhergehenden Detektor in den Flussweg anschließen.

#### Vorbereitungen

Siehe Flussverbindungen an der Messzelle ( Seite 71).

#### Gehen Sie wie folgt vor

Verbinden Sie die Einlasskapillare mit dem Einlass der Messzelle des Fluoreszenz-Detektors, wie es entsprechend des Systemaufbaus vorgegeben ist:

- Schließen Sie die Einlasskapillare direkt vom Säulenthermostaten an (siehe unten) –oder–
- Schließen Sie die Einlasskapillare vom Ausgang des UV-/VIS-Detektors an.
#### Verbinden der Einlasskapillare vom Säulenthermostaten

Verbinden Sie die Einlasskapillare zwischen dem Säulenthermostaten und dem Messzelleneinlass des Fluoreszenz-Detektors. Die Abbildung zeigt beispielhaft, wie die Verbindung vom Säulenthermostaten hergestellt werden kann.





1. Führen Sie die Einlasskapillare vom Säulenthermostaten durch das Führungsloch im Detektorgehäuse. Verwenden Sie das Führungsloch nahe dem Säulenthermostat.

**TIPP** Achten Sie stets darauf, die Kapillarverbindung zwischen Säulenthermostat und Messzelle möglichst kurz zu halten, um die Peak-Dispersion (d.h. Bandenverbreiterungseffekte durch Extra-Dispersionsvolumen) zu minimieren.

 Wenn sich der Säulenthermostat rechts vom Detektor befindet Führen Sie die Kapillare durch die untere Aussparung in der Zwischenwand. Stellen Sie sicher, dass Sie den geeigneten Zwischenwand-Einsatz für die Kapillare verwenden.

**TIPP** Kapillaren mit einem kleinen Außendurchmesser sind mit dem Einsatz mit Schlitz korrekt gesichert. Verwenden Sie den Dreheinsatz für Kapillaren, die nicht in den Einsatz mit Schlitz passen, wie isolierte Einlasskapillaren.

 Schließen Sie die Einlasskapillare an den Eingang der Messzelle an. Stellen Sie sicher, dass die Einlasskapillare angemessen in den Kapillarclip an der Vorderseite des Detektors gedrückt ist.

#### Verbinden der Einlasskapillare vom UV/VIS-Detektor

Schließen Sie die Kapillare zwischen dem Messzellenauslass des UV-/VIS-Detektors und dem Messzelleneinlass des Fluoreszenz-Detektors an. Die Abbildung zeigt beispielhaft, wie die Verbindung vom Diodenarray-Detektor hergestellt werden kann. Der Fluoreszenz-Detektor muss als letzter Detektor im Flussweg angeschlossen sein.

#### ACHTUNG

Beachten Sie die Rückdruckgrenze der Messzelle im Vanquish UV-/VIS-Detektor, welcher im Flussweg vor dem Fluoreszenz-Detektor verbunden ist. Schließen Sie die Kapillare vom UV-/VIS-Detektor direkt an den Einlass der Messzelle des Fluoreszenz-Detektors an. Schließen Sie keine zusätzlichen Komponenten im Flussweg zwischen den beiden Detektoren an. Informieren Sie sich in der *Betriebsanleitung für den UV-/ VIS-Detektor*.



Abbildung 24: Anschließen der Einlasskapillare vom Diodenarray-Detektor (Beispiel)

- Verbinden Sie die Einlasskapillare zwischen dem Säulenthermostaten und dem Messzelleneinlass des ersten Detektors. Folgen Sie den Anweisungen in der *Betriebsanleitung* für den ersten Detektor im Vanquish-System-Flussweg.
- 2. Führen Sie die Verbindungskapillare vom Auslass der Messzelle des ersten Detektors nach oben zum Fluoreszenz-Detektor.
- 3. Ziehen Sie den Kapillarclip an der Leckage-Wanne des Fluoreszenz-Detektors vorsichtig nach rechts, um den Clip zu öffnen.
- 4. Positionieren Sie die Kapillare hinter dem Kapillarclip.

5. Lassen Sie den Clip vorsichtig los und achten Sie darauf, die Kapillare nicht einzuklemmen. Die Kapillare muss sicher hinter dem Clip positioniert sein.



Abbildung 25: Fixieren der Kapillare hinter dem Kapillarclip an der Leckage-Wanne

Nr.	Beschreibung	Nr.	Beschreibung
1	Kapillare von der Messzelle	2	Kapillarclip

6. Schließen Sie die Einlasskapillare an den Messzellen-Einlass des Fluoreszenz-Detektors an. Stellen Sie sicher, dass die Einlasskapillare angemessen in den Kapillarclip an der Vorderseite des Detektors gedrückt ist.

#### 5.6.5.2 Anschließen der Detektor-Waste-Leitung

Vorbereitungen

1. Siehe Flussverbindungen an der Messzelle ( Seite 71).

#### Gehen Sie wie folgt vor

Schließen Sie die Detektor-Waste-Leitung zwischen dem Messzellenauslass des Fluoreszenz-Detektors und dem Abfall an. Die Abbildung zeigt die Schritte, um die Verbindung zum Messzellenausgang des Fluoreszenz-Detektors herzustellen.



Abbildung 26: Anschließen der Detektor-Waste-Leitung

- Schließen Sie die Waste-Leitung an den Messzellen-Auslass (OUT) (1) an.
- 2. Führen Sie die Waste-Leitung durch die obere Aussparung in der Zwischenwand (2).
- Führen Sie die Waste-Leitung durch die Schlauchführungen der Systemmodule unterhalb des Detektors zum Vanquish-Systemsockel (3).
- 4. Führen Sie die Waste-Leitung wie in der *Vanquish-System-Betriebsanleitung* beschrieben durch den dafür vorgesehenen Detektor-Waste-Auslass-Port im Systemsockel und anschließend vom Systemsockel zum Abfall.

#### TIPP

Die Waste-Leitung sollte direkt durch den Systemsockel zum Abfallbehälter gehen. Stellen Sie sicher, dass die Leitung gerade in den Schlauchführungen positioniert ist.

- 5. Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite des Schlauchs und dass die Schnitte nicht gequetscht sind.
- 6. Prüfen Sie die Waste-Leitung über den gesamten Flussweg: Stellen Sie sicher, dass die Waste-Leitung an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht ist.

#### 5.6.6 Ableiten von Flüssigkeiten in den Abfall

Flüssigkeiten aus Undichtigkeiten am Gerät sammeln sich in der Leckage-Wanne, wo sie durch den Drainage-Kanal rechts an der Leckage-Wanne in das Drainage-System geleitet werden.

Informationen dazu, wie Flüssigkeiten über die Drainage des Vanquish-Systems in den Abfall geleitet werden, finden Sie in der *Vanquish-System-Betriebsanleitung*.

### 5.7 Messen des Drucks in der Messzelle

Alle Module, Kapillaren und Waste-Leitungen nach der Messzelle tragen zum Druck in der Messzelle bei. Wenn Sie zusätzliche Module im Flussweg nach der Messzelle anschließen möchten, zum Beispiel Detektoren, Fraktionssammler oder Massenspektrometer, achten Sie darauf, dass der Druck in der Messzelle deren Druckspezifikation nicht übersteigt.

Wann

- Wenn Sie weitere Geräte und Kapillare im Flussweg nach der Messzelle anschließen möchten:
  - Bevor Sie die genannten zusätzlichen Geräte anschließen
  - Nachdem Sie die genannten weiteren Module angeschlossen haben
- Bevor Sie eine neue Methode verwenden
- Nachdem Sie Kapillaren im Flussweg nach der Messzelle ausgetauscht haben
- Abhängig von der Anwendung kann diese Prozedur bei der Fehlersuche helfen, zum Beispiel wenn Sie hohe Salzkonzentrationen verwenden.

#### Erforderliche Teile

- Verbindungsstück (zum Beispiel das Viper-Verbindungsstück aus dem System-Zubehör)
- Waste-Leitung und Kapillaren aus dem Detektor-Zubehör
- Eine der folgenden Lösungsmittelgemische mit hoher Viskosität:

Verwendung eines Gradienten	Lösungsmittelzusammensetzung
Wenn Sie Gradienten verwenden	Lösungsmittel für Ihre Anwendung. Verwenden Sie das Mischverhältnis mit der höchsten Viskosität.
Wenn Sie keine Gradienten verwenden	<ul> <li>60% Wasser und 40% Methanol</li> <li>–oder–</li> <li>70% Wasser und 30% Acetonitril</li> </ul>

#### Allgemeiner Ablauf der Prozedur

- Messen Sie den Rückdruck der Waste-Leitung (siehe Messen des Rückdrucks der Waste-Leitung (▶ Seite 80)). Flusspfad: Pumpe - Waste-Leitung
- Messen Sie den Rückdruck des Vanquish-Systems inklusive der Säule, der Detektor-Einlasskapillare und der Waste-Leitung (Messzelle ist nicht installiert) (siehe Messen des Rückdrucks des Vanquish-Systems (ohne Messzelle) (▶ Seite 81)). Flusspfad: Pumpe - Autosampler - Säule - Waste-Leitung
- Bestimmen Sie den Rückdruck der Transferkapillare und des zusätzlichen Moduls (siehe Bestimmen Sie den Rückdruck der Transferkapillare und des zusätzlichen Moduls (ohne Messzelle) (▶ Seite 83)).

Flusspfad: Pumpe - Autosampler - Säule - zusätzliches Modul (-Waste-Leitung)

- Bestimmen Sie den Rückdruck der Messzelle (siehe Bestimmen des Rückdrucks der Messzelle (▶ Seite 85)).
   Flusspfad: Pumpe - Autosampler - Säule - Messzelle - Waste-Leitung
- Berechnen Sie den Druck am Einlass der Messzelle (siehe Berechnen des Drucks am Einlass der Messzelle (▶ Seite 86)).
   Flusspfad: Pumpe - Autosampler - Säule - Messzelle - zusätzliches Modul (- Waste-Leitung)



#### WARNUNG—Austreten gefährlicher Substanzen aus Flussverbindungen

Fluss- und Kapillarverbindungen können mit Substanzen gefüllt sein, die ein Gesundheitsrisiko darstellen können. Lösungsmittel können herausspritzen, wenn Kapillaren bersten, aus ihren Fittingen rutschen, oder nicht korrekt festgezogen sind, oder wenn Kapillarverbindungen aus anderen Gründen offen sind.

- Tragen Sie eine geeignete Schutzausrüstung und folgen Sie der Guten Laborpraxis.
- Spülen Sie schädliche Substanzen vor Beginn der Wartungs- oder Servicearbeiten mit einem geeigneten Lösungsmittel aus.

#### Sehen Sie dazu auch

Messzellen (> Seite 178)

#### 5.7.1 Messen des Rückdrucks der Waste-Leitung

Im Folgenden wird beschrieben, wie der Rückdruck der Waste-Leitung (p0) gemessen wird.



Abbildung 27: Messen des Rückdrucks der Waste-Leitung

Nr.	Beschreibung	Nr.	Beschreibung
1	Detektor	3	Säule
2	Messzelle	4	Abfallbehälter
2a	Auslass der Messzelle	5	Zweiter Detektor
2b	Einlass der Messzelle	6	Waste-Leitung

In der folgenden Beschreibung wird davon ausgegangen, dass das Vanquish-System wie in der *System Betriebsanleitung* beschrieben aufgebaut wurde.

- 1. Schalten Sie das System ein (siehe Vanquish-System-Betriebsanleitung).
- 2. Lösen Sie die Einlass-Kapillare des Autosamplers vom Auslass der Pumpe/des statischen Mischers, falls zutreffend.
- 3. Schließen Sie die Waste-Leitung an den Auslass der Pumpe/des statischen Mischers an.
- 4. Führen Sie das andere Ende der Waste-Leitung in den Abfallbehälter.
- 5. Starten Sie den Pumpenfluss bei der maximalen Flussrate Ihrer Anwendung.Starten Sie den Pumpenfluss bei der maximalen Flussrate Ihrer Anwendung.

- 6. Prüfen Sie alle Flussverbindungen auf Undichtigkeiten:
  - Wenn eine Undichtigkeit auftritt: siehe Beheben von Undichtigkeiten (> Seite 171).
  - Wenn keine Undichtigkeit aufgetreten ist und sich der Druck stabilisiert hat: Fahren Sie mit dem nächsten Schritt fort.
- Wenn sich der Systemdruck stabilisiert hat, lesen Sie den Systemdruck in Chromeleon ab und notieren Sie sich den Wert für p0.

p0: Druckabfall der Waste-Leitung

- 8. Lösen Sie die Waste-Leitung vom Auslass der Pumpe/des statischen Mischers.
- 9. Schließen Sie die Einlass-Kapillare des Autosamplers am Auslass der Pumpe/des statischen Mischers an.

#### 5.7.2 Messen des Rückdrucks des Vanquish-Systems (ohne Messzelle)

Im Folgenden ist beschrieben, wie der Rückdruck des Vanquish-Systems inklusive der Säule, der Detektor-Einlasskapillare und Waste-Leitung (p1) gemessen wird.





Nr.	Beschreibung	Nr.	Beschreibung
1	Detektor	4	Verbindungsstück
2	Messzelle	5	Abfallbehälter
2a	Auslass der Messzelle	6	Zweiter Detektor
2b	Einlass der Messzelle	7	Waste-Leitung
3	Säule	8	Detektor-Einlasskapillare

- 1. Entfernen Sie die Detektor-Einlasskapillare vom IN-Port der Messzelle, falls zutreffend.
- 2. Entfernen Sie die Waste-Leitung vom OUT-Port der Messzelle, falls zutreffend.
- 3. Schließen Sie das freie Ende der Detektor-Einlasskapillare an das Verbindungsstück an.
- 4. Schließen Sie das freie Ende der Auslasskapillare an das andere Ende des Verbindungsstücks an.
- 5. Starten Sie den Pumpenfluss bei der maximalen Flussrate Ihrer Anwendung.
- 6. Prüfen Sie alle Flussverbindungen auf Undichtigkeiten:
  - Wenn eine Undichtigkeit auftritt: siehe Beheben von Undichtigkeiten (> Seite 171).
  - Wenn keine Undichtigkeit aufgetreten ist und sich der Druck stabilisiert hat: Fahren Sie mit dem nächsten Schritt fort.
- Wenn sich der Systemdruck stabilisiert hat, lesen Sie den Systemdruck in Chromeleon ab, und notieren Sie sich den Wert für p1.

p1: Druckabfall an der Säule, der Detektor-Einlasskapillare und der Waste-Leitung

8. Schalten Sie den Pumpenfluss ab.

#### Sehen Sie dazu auch

Flussverbindungen an der Messzelle (> Seite 71)

## 5.7.3 Bestimmen Sie den Rückdruck der Transferkapillare und des zusätzlichen Moduls (ohne Messzelle)

Im Folgenden ist zuerst beschrieben, wie der Rückdruck des zusätzlichen Moduls einschließlich der Säule, der Detektor-Einlasskapillare (und der Waste-Leitung) (p2) gemessen wird. Anschließend wird der Rückdruck der Transferkapillare und des zusätzlichen Moduls (p3) berechnet, indem der Rückdruck der Säule, der Detektor-Einlasskapillare (und der Waste-Leitung) (p1) abgezogen wird.



#### Abbildung 29: Messen des Rückdrucks eines zweiten Detektors (Beispiel)

Nr.	Beschreibung	Nr.	Beschreibung
1	Detektor	6	Auslass des zweiten Detektors
2	Messzelle	7	Einlass des zweiten Detektors
2a	Auslass der Messzelle	8	Zweiter Detektor
2b	Einlass der Messzelle	9	Waste-Leitung
3	Säule	10	Verbindungskapillare
4	Verbindungsstück	11	Detektor-Einlasskapillare
5	Abfallbehälter		

- 1. Schalten Sie das/die zusätzliche(n) Modul(e) im Flussweg nach der Messzelle ein. Informieren Sie sich in den *Betriebsanleitungen* der anderen Module.
- 2. Trennen Sie die Waste-Leitung vom Verbindungsstück.
- 3. *Nur wenn ein zweiter Detektor verwendet wird*: Schließen Sie die Waste-Leitung an den Auslass des zweiten Detektors an.

- 4. Verbinden Sie den Einlass des zusätzlichen Moduls mit dem freien Ende des Verbindungsstücks mit Hilfe der Verbindungskapillare, die bei der beabsichtigten Konfiguration für den Anschluss des zusätzlichen Moduls verwendet wird.
- 5. Starten Sie den Pumpenfluss bei der maximalen Flussrate Ihrer Anwendung.
- 6. Prüfen Sie alle Flussverbindungen auf Undichtigkeiten:
  - Wenn eine Undichtigkeit auftritt: siehe Beheben von Undichtigkeiten (> Seite 171).
  - Wenn keine Undichtigkeit aufgetreten ist und sich der Druck stabilisiert hat: Fahren Sie mit dem nächsten Schritt fort.
- Wenn sich der Systemdruck stabilisiert hat, lesen Sie den Systemdruck in Chromeleon ab, und notieren Sie sich den Wert für p2.

p2: Druckabfall an der Transferkapillare und am zusätzlichen Modul einschließlich Säule, Detektor-Einlasskapillare und Waste-Leitung

- 8. Schalten Sie den Pumpenfluss ab.
- 9. Berechnen Sie die Differenz zwischen den zwei gemessenen Druckwerten: p3 = p2 - p1.
  p3: Druckabfall an der Transferkapillare und am zusätzlichen Modul

p2: Druckabfall an der Transferkapillare und am zusätzlichen Modul einschließlich Säule, Detektor-Einlasskapillare und Waste-Leitung

p1: Druckabfall an der Säule, der Detektor-Einlasskapillare und der Waste-Leitung

10. Vergleichen Sie p3 mit dem Druck-Grenzwert der Messzelle.

Situation	Schritte
Wenn p3 deutlich unterhalb des Druck- Grenzwerts der Messzelle liegt	Fahren Sie mit Bestimmen des Rückdrucks der Messzelle (▶ Seite 85) fort.
Wenn p3 nahe am Druck- Grenzwert für die Messzelle liegt oder darüber	Verringern Sie schrittweise die Flussrate und wiederholen Sie die Messung bis p3 deutlich unterhalb des Druck-Grenzwerts der Messzelle liegt. –oder–
	Verbindungskapillare) und führen Sie die Prozedur erneut aus, bis p3 deutlich unterhalb des Druck- Grenzwerts der Messzelle liegt.

#### Sehen Sie dazu auch

Messzellen (> Seite 178)

#### 5.7.4 Bestimmen des Rückdrucks der Messzelle

Im Folgenden wird zuerst beschrieben, wie der Rückdruck der Messzelle (p4) einschließlich der Säule, der Detektor-Einlasskapillare und der Waste-Leitung gemessen wird. Anschließend wird der Rückdruck der Säule, der Detektor-Einlasskapillare und der Waste-Leitung (p1) abgezogen.



Abbildung 30: Messen des Rückdrucks der Messzelle einschließlich der Säule, der Detektor-Einlasskapillare und der Waste-Leitung

Nr.	Beschreibung	Nr.	Beschreibung
1	Detektor	3	Detektor-Einlasskapillare
2	Messzelle	4	Abfallbehälter
2a	Auslass der Messzelle	5	Waste-Leitung
2b	xEinlass der Messzelle		

- 1. Stellen Sie die Flussverbindungen wie unter Flussverbindungen an der Messzelle (▶ Seite 71) beschrieben her.
- 2. Starten Sie den Pumpenfluss bei der Flussrate Ihrer Anwendung.
- 3. Prüfen Sie alle Flussverbindungen auf Undichtigkeiten:
  - Wenn eine Undichtigkeit auftritt: siehe Beheben von Undichtigkeiten (> Seite 171).
  - Wenn keine Undichtigkeit aufgetreten ist und sich der Druck stabilisiert hat: Fahren Sie mit dem nächsten Schritt fort.
- Wenn sich der Systemdruck stabilisiert hat, lesen Sie den Systemdruck in Chromeleon ab, und notieren Sie sich den Wert für p4.

p4: Druckabfall an der Messzelle einschließlich Säule, Detektor-Einlasskapillare und Waste-Leitung

5. Schalten Sie den Pumpenfluss ab.

6. Berechnen Sie die Differenz zwischen den zwei gemessenen Druckwerten: p5 = p4 - p1.
p5: Druckabfall an der Messzelle

p4: Druckabfall an der Messzelle einschließlich Säule, Detektor-Einlasskapillare und Waste-Leitung

p1: Druckabfall an der Säule, der Detektor-Einlasskapillare und der Waste-Leitung

#### 5.7.5 Berechnen des Drucks am Einlass der Messzelle

Im Folgenden wird beschrieben, wie der Rückdruck am Einlass der Messzelle (p6) für die gewünschte Konfiguration ermittelt wird.



Abbildung 31: Errechnen des Rückdrucks am Messzelleneinlass in der angestrebten Systemkonfiguration (hier mit zweitem Detektor)

Nr.	Beschreibung	Nr.	Beschreibung
1	Detektor	5	Abfallbehälter
2	Messzelle	6	Auslass des zweiten Detektors
2a	Auslass der Messzelle	7	Einlass des zweiten Detektors
2b	Einlass der Messzelle	8	Zweiter Detektor
3	Detektor-Einlasskapillare	9	Waste-Leitung
4	Säule	10	Verbindungskapillare

- Berechnen Sie P6 = P5 + P3 + p0. p6: Druck am Messzelleneingang
  - p5: Druckabfall an der Messzelle
  - p3: Druckabfall an der Transferkapillare und am zusätzlichen Modul
  - p0: Druckabfall an der Waste-Leitung

2. Vergleichen Sie p6 mit dem Druck-Grenzwert der Messzelle.

Situation	Schritte
p6 liegt mindestens 5 bar unterhalb des Druck- Grenzwerts der Messzelle	Die Konfiguration kann verwendet werden.
p6 liegt weniger als 5 bar unterhalb des Druck- Grenzwerts der Messzelle oder liegt über dem Druck-Grenzwert	Betreiben Sie die Systemkonfiguration bei einer niedrigeren Flussrate. –oder– Verändern Sie die Systemkonfiguration (zum Beispiel die Verbindungskapillare), damit die Druckspezifikation des Messzelle erreicht werden kann.

#### Sehen Sie dazu auch

Messzellen (> Seite 178)

### 5.8 Einschalten des Gerätes

#### TIPP

Bevor Sie ein Modul des Vanquish-Systems zum ersten Mal einschalten vergewissern Sie sich, dass die Chromatographie-Software auf dem Datensystemrechner installiert ist. Die erforderlichen USB-Treiber werden automatisch geladen und das Windows™-Betriebssystem kann das Gerät erkennen, wenn dieses eingeschaltet ist.

Schalten Sie das Gerät wie folgt ein:

- Vergewissern Sie sich, dass der Ein-/Ausschalter vorne links am Vanquish-Systemsockel (System-Ein-/Ausschalter) gedrückt ist. Wenn der Ein-/Ausschalter heraussteht, drücken Sie den Ein-/ Ausschalter, um den Systemsockel einzuschalten.
- 2. Schalten Sie das Gerät über den Hauptnetzschalter ein.

Schalten Sie das Gerät über den Hauptnetzschalter aus, wenn Sie dazu aufgefordert werden, zum Beispiel für Wartungsarbeiten. Das Drücken des System-Ein-/Ausschalters reicht nicht aus, um das Gerät vollständig auszuschalten.

#### Sehen Sie dazu auch

Ein- und Ausschalten (▶ Seite 97)

### 5.9 Einrichten des Gerätes in der Software

Diese Anleitung setzt voraus, dass die Chromatographie-Software bereits auf dem Datensystemrechner installiert und eine gültige Lizenz verfügbar ist.

Weitere Informationen zum Einrichten des Vanquish-Systems in der Software, finden Sie in der *Vanquish-System-Betriebsanleitung*.

Genauere Informationen zu den Einstellungen auf den einzelnen Konfigurationsseiten erhalten Sie in der Hilfe der verwendeten Software.

# **6 Betrieb**

Dieses Kapitel enthält Informationen zu den Bedienelementen des Gerätes, zum Routinebetrieb und zur Außerbetriebnahme.

### 6.1 Einführung in dieses Kapitel

Die Informationen in diesem Kapitel setzen voraus, dass die Ersteinrichtung des Gerätes bereits abgeschlossen ist. Ist dies nicht der Fall, folgen Sie den Anweisungen in Installation (▶ Seite 47) bevor Sie fortfahren.

Eine kurze Beschreibung der Gerätesteuerung und automatisierten Probenanalyse mit der Software Chromeleon finden Sie in der Vanquish-System-Betriebsanleitung. Einzelheiten zu Steuerbefehlen und Bedienung des Geräts finden Sie in der Chromeleon-Hilfe.

Die Software-Beschreibungen in diesem Anleitung beziehen sich auf Chromeleon 7. Die Terminologie kann sich gegebenenfalls geringfügig von der Terminologie anderer Software-Versionen unterscheiden.

### 6.2 Sicherheitshinweise zum Betrieb

Beachten Sie beim Betrieb des Gerätes folgende Sicherheitshinweise:



Beachten Sie alle Warnhinweise und Vorsichtsmaßnahmen in Sicherheitsmaßnahmen (▶ Seite 21).



#### VORSICHT—Heiße Oberflächen

Oberflächen im Inneren des Geräts können sich während des Betriebs erhitzen. Das Berühren von heißen Komponenten kann zu Verbrennungen führen.

Warten Sie, bis die heißen Oberflächen abgekühlt sind, bevor Sie diese berühren.

#### ACHTUNG

Beachten Sie auch folgende Hinweise:

Wenn der Pumpenfluss unterbrochen wird, ergreifen Sie geeignete Maßnahmen, um die Messzelle zu schützen. Beachten Sie die Hinweise zum Gebrauch von Messzellen in Hinweise zum Gebrauch von Messzellen (> Seite 100).

### 6.3 Bedienelemente

Das Gerät wird hauptsächlich über einen Rechner gesteuert, auf dem die Chromatographie-Software installiert ist.

Zusätzlich stehen folgende Bedienelemente am Gerät zur Verfügung:

- Tastatur Über die Tasten können Sie einige Funktionen direkt am Gerät ausführen.
- Statusanzeigen Die LEDs (Light Emitting Diodes) der LED-Leiste (Statusanzeige) an der Vorderseite des Geräts sowie die **STATUS**-LED auf der Tastatur ermöglichen einen schnellen, visuellen Überblick über den Betriebszustand des Geräts.

#### 6.3.1 Tastatur

Über die Tastatur im Gerät können Sie einige Funktionen direkt am Gerät ausführen. Beim Drücken eines Knopfes bestätigt ein akustisches Signal, dass die Aktion ausgeführt wird.



Abbildung 32: Tastatur

#### STATUS

Über die **STATUS** LED können Sie schnell den Betriebszustand des Gerätes erkennen.

Wenn die Gerätetüren geschlossen sind, zeigt die LED-Leiste an der Vorderseite den Betriebszustand an.

Weitere Informationen zur Statusanzeige finden Sie in Statusanzeigen (> Seite 95).

#### MUTE ALARM

Ein akustisches Signal ertönt, wenn das Gerät ein Problem erkannt hat, zum Beispiel eine Undichtigkeit. Drücken Sie diese Taste, um das akustische Signal für den aktuellen Alarm auszuschalten. Beseitigen Sie die Ursache für die Undichtigkeit innerhalb von 10 Minuten. Andernfalls ertönt das Signal erneut. Wenn das Gerät ein anderes Problem erkennt, ertönt sofort wieder ein akustisches Signal.

#### INIT

Durch Drücken der INIT-Taste können Sie eine Initialisierung des Geräts durchführen, wobei die Startpositionen der Gittermotoren und des Filterrads bestimmt werden. Die LED neben der Taste zeigt den Initialisierungsstatus an:

LED	Beschreibung
Aus (dunkel)	Der Detektor ist nicht initialisiert, oder ausgeschaltet.
Grün, blinkend	Der Detektor initialisiert gerade.
Grün	Der Detektor ist initialisiert.

Wenn die LED nach der Initialisierung erlischt, sollten Sie die korrekte Installation der Messzelle überprüfen und anschließend die **INIT**-Taste erneut drücken.

#### 6.3.2 Statusanzeigen

Die LED-Statusleiste an der Vorderseite des Geräts und die **STATUS**-LED an der Tastatur im Innenraum informieren Sie über den Zustand des Geräts.

#### LED-Leiste

Die LED-Leiste zeigt die Informationen an, wenn das Gerät geschlossen ist. Ist das Gerät in der Software Chromeleon verbunden, zeigt die LED-Leiste gegebenenfalls weniger Informationen an.

LED-Leiste	Beschreibung
Aus (dunkel)	Das Gerät ist ausgeschaltet.
Gedimmt	Die Gerätetüren sind geöffnet.
Gelb, langsam blinkend	Das Gerät ist eingeschaltet, jedoch nicht in der Software Chromeleon verbunden.
Gelb	Das Gerät ist in der Software Chromeleon verbunden, es ist aber nicht initialisiert.
Grün, blinkend	Das Gerät initialisiert gerade.
Grün	Das Gerät ist initialisiert, es läuft jedoch keine Datenaufnahme.
Blau, Lauflicht	Eine Datenaufnahme läuft.
Blau	Eine Injektion oder Sequenz läuft.
Rot	Ein Problem oder Fehler ist aufgetreten. Prüfen Sie den Chromeleon Audit Trail auf eine entsprechende Meldung. Abhilfemaßnahmen finden Sie unter Fehlersuche (> Seite 163).

#### STATUS LED

### Die **STATUS** -LED auf der Tastatur im Gerät zeigt die folgenden Informationen an:

STATUS LED	Beschreibung
Aus (dunkel)	Das Gerät ist ausgeschaltet.
Grün	Das Gerät funktioniert ordnungsgemäß.
Rot	Ein Problem oder Fehler ist aufgetreten. Prüfen Sie den Chromeleon Audit Trail auf eine entsprechende Meldung. Abhilfemaßnahmen finden Sie unter Fehlersuche (> Seite 163).

### 6.4 Ein- und Ausschalten

Der Netzschalter am Gerät ist der Hauptnetzschalter zum Ein- und Ausschalten des Gerätes. Der Hauptnetzschalter wird bei der Inbetriebnahme des Gerätes eingeschaltet.

Für eine einfachere Bedienung können Sie den Schalter vorne links am Vanquish-Systemsockel (System-Ein-/Ausschalter) zum Ein- und Ausschalten verwenden.

Beachten Sie Folgendes:

- *Alle* Module des Vanquish-Systems, die über einen System-Interlink-Port mit dem Systemsockel verbunden sind, werden durch Drücken des System-Ein/Ausschalters gleichzeitig ein- oder ausgeschaltet.
- Der Sockel ist eingeschaltet, wenn der System-Ein-/Ausschalter gedrückt ist. Der Sockel ist ausgeschaltet, wenn der System-Ein-/ Ausschalter heraussteht.
- Wenn der Hauptnetzschalter eines Geräts ausgeschaltet ist, können Sie das Gerät nicht über den System-Ein- /Ausschalter einschalten.
- Um ein Gerät vollständig auszuschalten, müssen Sie es über den Hauptnetzschalter des Geräts ausschalten. Das Drücken des System-Ein-/Ausschalters reicht nicht aus, um das Gerät vollständig auszuschalten.

Beim Einschalten des Gerätes ist Folgendes zu beobachten, wenn die Gerätetüren geschlossen sind:

- Das Gerät führt einen Selbsttest durch. (Der Test dauert circa 30 Sekunden.) Während des Selbsttests werden alle wesentlichen Baugruppen auf korrekte Funktion überprüft.
- Nach dem Selbsttest beginnt der Initialisierungsvorgang. Während dieser Zeit blinkt die LED-Leiste grün.
- Nach erfolgreicher Initialisierung leuchtet die LED-Leiste grün.
   Wenn die LED-Leiste rot leuchtet, sollten Sie überprüfen, ob die Messzelle korrekt eingebaut ist und anschließend die INIT-Taste auf der Tastatur drücken, um eine erneute Initialisierung zu starten. Die LED-Leiste und die LED neben der INIT-Taste leuchten nach erfolgreicher Initialisierung grün.

### 6.5 Vorbereiten des Gerätes für den Betrieb

In diesem Abschnitt finden Sie weitere Schritte, die zur Vorbereitung des Gerätes für den Betrieb und die Probenanalyse erforderlich sind.

Vor der ersten Inbetriebnahme des Gerätes

Bereiten Sie das Gerät für die Erstinbetriebnahme vor; beachten Sie dabei Folgendes:

#### ACHTUNG

Spülen Sie den System-Flussweg gründlich, bevor Sie das Gerät zum ersten Mal in Betrieb nehmen:

- Wenn Sie Geräte oder Komponenten im System installieren, spülen Sie diese immer in den Abfall, bevor Sie diese in den Flussweg des Systems aufnehmen. Folgen Sie den Anweisungen zum Spülen der Vanquish-Module in der Vanquish-System-Betriebsanleitung.
- Wenn Sie das Gerät mit einer Messzelle betreiben, die zuvor gelagert wurde, könnte die Messzelle mit Lösungsmittel gefüllt sein. Verwenden Sie entsprechend Lösungsmittel, die mit diesem Lösungsmittel mischbar sind, oder verwenden Sie einen geeigneten Löslichkeitsvermittler.

#### Vor dem Beginn einer Probenanalyse

Bevor Sie mit einer Probenanalyse beginnen:

- Überprüfen Sie den Füllstand in den Lösungsmittelbehältern.
   Vergewissern Sie sich, dass die Lösungsmittelmenge für die Analyse ausreicht.
- Vergewissern Sie sich, dass die Gerätetüren aller Module im Vanquish-System geschlossen sind.
- Beachten Sie die Hinweise zum Gebrauch von Messzellen in Hinweise zum Gebrauch von Messzellen (▶ Seite 100).
- Stellen Sie sicher, dass das Chromatographie-System ausreichend äquilibriert ist (siehe weiter unten).

#### System-Äquilibrierung

Die System-Äquilibrierung sollte Folgendes umfassen:

• Spülen *aller* Kanäle der Pumpe (auch der Kanäle, die für die Anwendung nicht verwendet werden)

- Spülen des gesamten Chromatographie-Systems mit dem Anfangseluenten, um Lösungsmittel der vorherigen Analyse auszuspülen
- Heizen (oder Kühlen) aller temperaturgesteuerten Geräte im System auf die Anfangstemperatur. Temperaturgesteuerte Geräte können zum Beispiel sein
  - Säulenthermostat und Nachsäulenwärmetauscher
  - Thermostatisierter Probenraum im Autosampler
  - Messzelle in einem Fluoreszenz-Detektor
  - Verdampfungsrohr in einem Charged-Aerosol-Detektor
- Einschalten der Lampe (oder Lampen) im UV-/VIS-Detektor
- Beobachten des Pumpendrucks und der Druckpulsation sowie überprüfen, dass der Druck stabil ist und die Pulsation in einem vernünftigen Bereich für die Anwendung liegt
- Beobachten des Detektorsignals und überprüfen, ob das Detektorsignal stabil ist, so dass Drift und Signalrauschen in einem vernünftigen Bereich für die Anwendung liegen
- Durchführen eines Autozero der Detektor-Basislinie

**TIPP** Die Software Chromeleon unterstützt Prozeduren, um ein Chromatographie-System automatisch in der Software zu starten (**Smart Startup**). Der Startvorgang beinhaltet Prozeduren für die System-Äquilibrierung. Einzelheiten hierzu finden Sie in der *Chromeleon-Hilfe*.

### 6.6 Hinweise zum Gebrauch von Messzellen

#### ACHTUNG

Die Messzellen reagieren empfindlich auf Beschädigung, Verschmutzung und Staub.

- Behandeln Sie Messzellen mit Vorsicht. Siehe Hinweise zum Gebrauch von Messzellen (▶ Seite 141).
- Beachten Sie bei der Bedienung des Detektors die unten stehenden Hinweise.

#### Betriebsbedingungen

Beachten Sie den spezifizierten maximalen Druck-Grenzwert für die Messzelle. Siehe Spezifikationen für die Messzellen unter Messzellen (**b** Seite 178).

#### Fehlersuche bei den Messzellen

Tritt eine Undichtigkeit in einer Messzelle auf, stoppen Sie den Pumpenfluss, entfernen Sie die Messzelle möglichst schnell aus dem Detektor, entfernen Sie ausgetretene Flüssigkeiten im Messzellenschacht gründlich und tauschen Sie die Messzelle aus.

#### Betriebsunterbrechung

Wenn der Pumpenfluss unterbrochen wird, ergreifen Sie geeignete Maßnahmen, um die Messzelle zu schützen.

- Halten Sie die Datenaufnahme an, falls diese noch läuft.
- Lassen Sie keine Substanzen, insbesondere aggressive Lösungsmittel, für längere Zeit ohne Fluss in der Messzelle stehen.

**TIPP** Um bei Betriebsunterbrechungen Schäden an der Detektoroptik durch Staubpartikel zu vermeiden, achten Sie darauf, dass am Detektor immer eine Messzelle oder die mitgelieferte Messzellen-Abdeckung installiert ist.

#### Aufbewahrung

- Um Algenbildung zu vermeiden sollte die Messzelle nicht mit reinem Wasser gefüllt sein. Geben Sie zum Beispiel 10%iges Isopropanol in HPLC-Qualität zu.
- Verschließen Sie den Einlass und Auslass der Messzelle mit den Verschlussstopfen, die bei Auslieferung an der Messzelle installiert waren. Wenn Sie andere Stopfen verwenden und diese zudrehen, kann dies die Messzelle zerstören.

### 6.7 Betriebsmodi des Detektors

Die folgenden fünf Betriebsmodi stehen zur Verfügung:

- Single-Channel-Modus
- Multi-Channel-Modus
- Zero-Order-Modus
- Single Spectrum Scan (Erfassung von Einzelspektren)
- FL Field Acquisiton

#### 6.7.1 Single-Channel-Modus

Ein Fluoreszenz-Detektor wird üblicherweise im Single-Channel-Modus betrieben, das heißt, es wird *ein* Anregungs- / Emissionswellenlängenpaar über die Zeit gemessen. Andere Wellenlängenpaare und Detektionsparameter können jederzeit während der Probentrennung eingestellt werden, um sie an die Besonderheiten der Analyten anzupassen. Diese Anpassung sollte bevorzugt dann erfolgen, wenn kein Peak detektiert wird. Da hier die höchstmöglichen Datenraten und das beste Signal-Rausch-Verhältnis erreicht werden können, verwenden Sie bei sehr schmalen Peaks stets den Single-Channel-Modus.

#### 6.7.2 Multi-Channel-Modus (Mehrkanalbetrieb, nur VF-Detektoren)

Im Multi-Channel-Modus werden beide Monochromatoren schnell zwischen den ausgewählten Wellenlängen hin- und her bewegt. Es ist daher eine einzige Analyse ausreichend, um auf mehreren Kanälen zu messen. Der Detektor kann bis zu vier Wellenlängenpaare gleichzeitig messen. Beachten Sie jedoch Folgendes:

- Jeder zusätzliche Kanal erhöht die Zeit, die der Detektor zum Einstellen der Gitterpositionen benötigt, reduziert die maximal erreichbare Datenaufnahmerate und führt zu einem höheren Basislinienrauschen.
- Ein ständiges Umschalten zwischen den Wellenlängen führt über einen längeren Zeitraum zum Verschleiß. Der Multi-Channel-Modus wird daher nur für die Methodenentwicklung empfohlen, jedoch nicht für den dauerhaften Routinebetrieb.

 Alternativ können Sie die Wellenlängen nach Bedarf zwischen den einzelnen Peaks auf einem einzigen Datenkanal mit Hilfe der Software Chromeleon umschalten. Zusätzlich zur Anregungs- und Emissionswellenlänge können Sie die Empfindlichkeit, den Emissionsfilter, den verwendeten PMT sowie den Lampenmodus schalten.

#### Einstellen der Multi-Channel Performance (Messqualität)

Im Multi-Channel-Modus können Sie die Messqualität (Multi-Channel Performance) festlegen, welche bestimmt, ob eine Messung schneller, dafür aber mit mehr Rauschen durchgeführt wird, oder ob möglichst wenig Rauschen gewünscht wird, was eine längere Messzeit und daher geringere Datenaufnahmerate bedeutet. Mögliche Werte sind: **UltraFast, Fast, Standard, LowNoise, UltraLowNoise**.

Die Datenaufnahmerate und Ansprechzeit werden im Multi-Channel-Modus automatisch anhand der gewählten Multi-Channel Performance und der weiteren Einstellungen festgelegt und bei Wellenlängenschaltungen während der Messung entsprechend angepasst.

#### Aufzeichnen von mehreren Kanälen in der Software Chromeleon

Verwenden Sie den Instrument Method Wizard (Modus Advanced), um die Parametereinstellungen für alle Kanäle sowie die Multi-Channel Performance (Messqualität) festzulegen.

#### Sehen Sie dazu auch

Optimale Anregungs- und Emissionswellenlänge bestimmen
 () Seite 110)

#### 6.7.3 Zero-Order-Modus

Im Zero-Order-Modus reflektiert das Gitter des Emissions-Monochromators das gesamte Emissionsspektrum der Probe auf den PMT, nicht nur eine einzelne Wellenlänge. Der Anregungs-Monochromator wird wie gewohnt auf eine einzelne Wellenlänge eingestellt.

Verwenden Sie diesen Modus in der Methodenentwicklung, wenn Sie die Retentionszeiten und Emissionswellenlängen der verschiedenen Substanzen in Ihrer Probe nicht kennen. Da der gesamte Bereich der Emissionswellenlängen erfasst wird, können die Retentionszeiten aller Substanzen in einem einzigen Durchlauf erfasst werden, solange alle Substanzen bei der gewählten Anregungswellenlänge angeregt werden.

#### TIPP

Mit Hilfe der Filterrad-Einstellung (nur VF-Detektoren) können Sie unerwünschte Wellenlängen ausblenden. Stellen Sie die Filterrad-Einstellung zum Beispiel auf eine Wellenlänge von 370 nm, um den Emissionswellenlängenbereich auf Wellenlängen oberhalb der Filterwellenlänge zu beschränken.

Der Zero-Order-Modus eignet sich außerdem für Proben, die Licht mit einer ungewöhnlich großen Bandbreite emittieren. In diesem Fall ist die gemessene Intensität im Zero-Order-Modus höher als im Normalbetrieb, wo der Großteil des emittierten Lichts vom Emissions-Monochromator verworfen wird. Hiermit können Sie gegebenenfalls eine bessere Detektionsgrenze erreichen.

Ein Beispiel zur Methodenentwicklung mit dem Zero-Order-Modus finden Sie im Fluorescence Method Development Handbook.

#### TIPP

Setzen Sie zum Aktivieren des Zero-Order-Modus in der Software Chromeleon den Wert für **EmWavelength** auf **ZeroOrder**.

#### Sehen Sie dazu auch

Filterrad (Nur VF-Detektoren) ( Seite 115)

#### 6.7.4 Single Spectrum Scan (Erfassung von Einzelspektren)

Bei einem Single Spectrum Scan werden der Anregungs-Monochromator oder der Emissions-Monochromator (oder auch beide simultan) über einen einstellbaren Wellenlängenbereich durchgefahren, während die Intensität des Fluoreszenz-Signals in Abhängigkeit von der Wellenlänge kontinuierlich gemessen und aufgezeichnet wird. Nehmen Sie zum Beispiel Spektren auf, um die optimalen Emissions- und Anregungswellenlängen zu bestimmen. Es stehen drei Arten von Scanmodi zur Verfügung: Excitation Scan, Emission Scan und Synchronous Scan.

Um einen stabilen Zustand in der Messzelle des Fluoreszenz-Detektors zu erreichen sollte während der Scans der Pumpenfluss angehalten oder auf eine sehr geringe Flussrate eingestellt werden. Wenn der Scan während der Elution eines Peaks abgeschlossen werden kann und der Scan wesentlich schneller als die zeitliche Halbwertsbreite des Peaks ist, kann der Pumpenfluss konstant gehalten werden. Weitere Informationen finden Sie in der *Chromeleon-Hilfe*.

#### Aufzeichnen eines Basislinienspektrums

		Für alle Scan-Verfahren kann ein Basislinienspektrum mit den selben Parametern aufgenommen werden. Dieses wird in der Software Chromeleon gespeichert und automatisch von den anderen Spektren subtrahiert. Das Ergebnis ist ein Differenzspektrum, d.h. die Autofluoreszenz des verwendeten Lösungsmittels wird eliminiert. Beachten Sie, dass Chromeleon eine Warnung ausgibt, wenn das Basislinienspektrum fehlt.
		Werden die Parameter für einen Scan geändert, wird das Basislinienspektrum automatisch gelöscht und muss neu aufgenommen werden. Sie können zudem jederzeit ein bestehendes Basislinien- spektrum über den Befehl <b>ClearBaseLine</b> löschen.
	Excitation Scan	
		Die Wellenlänge auf dem Emissions-Monochromator wird konstant gehalten, während der Anregungs-Monochromator einen Wellenlängenbereich scannt. Sie erhalten das Anregungsspektrum der Probe.
	Emission Scan	
		Die Wellenlänge auf dem Anregungs-Monochromator wird konstant gehalten, während der Emissions-Monochromator einen Wellenlängenbereich scannt. Sie erhalten das Emissionsspektrum der Probe.
	Synchronous Scan	
		Ein festgelegter Bereich der Anregungswellenlänge wird gescannt, während synchron dazu mit einem benutzerdefinierten Offset die Emissionswellenlänge gescannt wird. Dies ermöglicht Ihnen, ein geeignetes Wellenlängenpaar für erste Experimente zu bestimmen. Die optimalen Anregungs- und Emissionswellenlängen können jedoch nur mit separaten Excitation- und Emission-Scans ermittelt werden.
So führen Sie einen Single Spectrum Scan durch		
		Informationen zu Spektrenscans und Beispiele für Instrumentenmethoden finden Sie in der <i>Chromeleon-Hilfe</i> .



#### VORSICHT

- Während des Scan-Vorgangs können keine anderen Befehle an den Fluoreszenz-Detektor gesendet werden.
- Die Dauer des Scan-Vorgangs hängt von der Größe des zu scannenden Bereiches sowie der gewählten Scan-Geschwindigkeit ab.
- Wählen Sie vor jedem Scan über den Parameter ScanSensitivity eine geeignete Detektorempfindlichkeit. Wenn die Detektorempfindlichkeit zu hoch gewählt ist, wird kein Spektrum gespeichert und eine Meldung erscheint im Audit Trail.
- In der Nähe von starken Spektraländerungen solche zeigen sich insbesondere, wenn nahe der Anregungswellenlänge gescannt wird – können beim Scannen Artefakte entstehen, zum Beispiel im Raman-Emissionsspektrum von Wasser nahe der Anregung von 350 nm. Wählen Sie in diesem Fall einen größeren Abstand des Wellenlängenbereichs von der Flanke (größer als der minimale Abstand von 20 nm). Wählen Sie für den Raman-Emissionsscan zum Beispiel eine Startwellenlänge von 385 nm.

#### 6.7.5 FL Field Acquisition

Mit Hilfe des Modus FL Field Acquisiton können Sie die Retentionszeiten und Emissionsmaxima bestimmen. Im Gegensatz zu einem Single Spectrum Scan wird der gewählte Spektrenbereich dabei nicht *einmal*, sondern kontinuierlich gescannt, so dass ein Spektrenfeld ähnlich dem 3D-Feld eines Diodenarray-Detektors entsteht. Folgende Modi stehen zur Verfügung:

- Excitation (Anregung)
- Emission
- Synchronous

Es ist nicht möglich, während der Aufzeichnung eines FL-Feldes im Modus FL Field Acquisiton ein Chromatogramm aufzunehmen.

Der Modus FL Field Acquisiton ist nur in der Software Chromeleon 7 möglich. Verwenden Sie ein 3D-Datenfeld, um die Retentionszeiten sowie die optimalen Emissions- oder Anregungswellenlängen zu bestimmen. Weitere Informationen zu 3D-Datenfeldern finden Sie in der *Chromeleon-Hilfe*.

Ein Beispiel zur Methodenentwicklung mit dem Modus FL Field Acquisiton finden Sie im Fluorescence Method Development Handbook.

### 6.8 Wichtige Einstellungen für den Betrieb

Die Kommandos und Parameter in dieser Tabelle sollten für den einfachen Routinebetrieb des Geräts berücksichtigt werden. Sie können diese Einstellungen in der Regel über die Benutzeroberfläche in Chromeleon öffnen.

Ist einer der unten genannten Parameter nicht in der Software Chromeleon vorhanden, aktualisieren Sie gegebenenfalls die Firmwareund Chromeleon-Version.

**TIPP** Der Instrument Method Wizard bietet verschiedene Parameter-Ansichten für den Detektor. Abhängig von den einzustellenden Parametern kann die gewünschte Ansicht (zum Beispiel **Easy** oder **Advanced**) ausgewählt werden.

Weitere Informationen finden Sie in der *Chromeleon-Hilfe und Benutzerdokumentation*.

Parameter	Beschreibung
Emissionsparameter	Die Emissions-Kanäle ( <b>Emission_x</b> ) sind die Signalkanäle für die Aufnahme eines Anregungs-/ Emissions- Wellenlängenpaares. Die Signalparameter sind für jeden Signalkanal separat einstellbar. Hinweise zur Auswahl der Signalparameter finden Sie unter Optimieren der Geräteleistung (▶ Seite 108).
Anregungs- wellenlänge	Legt die Anregungswellenlänge in nm fest. Wählen Sie die Anregungswellenlänge auf dem Absorptionsmaximum des zu analysierenden Probenbestandteils.
Emissions- wellenlänge	Legt die Emissionswellenlänge in nm fest. Die Emissionswellenlänge muss mindestens 20 nm über der Anregungswellenlänge liegen. <b>ZeroOrder</b> : Das Gitter des Emissions-Monochromators wird auf die Position der 0. Ordnung eingestellt. Einzelheiten finden Sie unter Zero-Order-Modus (▶ Seite 102).
Empfindlichkeit	Legt die Empfindlichkeit des Detektors fest, siehe Sensitivity (Detektorempfindlichkeit) (▶ Seite 111).
Datenaufnahmerate	Legt fest, wie viele Datenpunkte die Software Chromeleon pro Sekunde (Hz) vom Detektor aufnimmt und als Rohdaten abspeichert. Weitere Informationen finden Sie unter Datenaufnahmerate und Ansprechzeit (Single-Channel- Modus) (> Seite 117).
Lampen-Modus	Legt den Betriebsmodus für die Xenon-Blitzlampe fest. Eine höhere Frequenz verbessert das Basislinienrauschen und führt damit zu einer höheren Empfindlichkeit. Eine geringere Frequenz verlängert die Lebensdauer der Lampe. Einzelheiten finden Sie in Lampen-Modus (> Seite 119).

Parameter	Beschreibung
Betriebsdauer (Effective Lamp Age)	Liefert Informationen über die Betriebsdauer der Lampe; diese Funktion soll helfen, festzustellen, wann eine Lampe ausgetauscht werden muss. Weitere Informationen finden Sie unter Überwachen der Lampenbetriebsdauer () Seite 136).
Initialize	Führt eine Initialisierung des Geräts durch, wobei die Startpositionen der Gittermotoren und des Filterrads bestimmt werden. Führen Sie den Befehl aus, wenn beim Einschalten des Detektors die Meldung "Not initialized" im Chromeleon Audit Trail erscheint.
Leakerkennung	Bei Auslieferung des Geräts ist die Leakerkennung standardmäßig aktiviert ( <b>Leak Sensor Mode = Enabled</b> ). Dies ist die bevorzugte Einstellung.

### 6.9 Optimieren der Geräteleistung

In diesem Abschnitt finden Sie Informationen zur höchsten Leistungsfähigkeit des Gerätes und Hinweise, wie Sie die Leistung noch weiter optimieren können.

#### Fluorescence Method Development Handbook

Grundlegende Informationen zur Optimierung der Detektorleistung erhalten Sie in den folgenden Kapiteln dieser Betriebsanleitung. Das Fluorescence Method Development Handbook, das (in englischer Sprache) mit dem Gerät ausgeliefert wird, führt durch die verschiedenen Schritte der Methodenentwicklung und Optimierung, gibt Empfehlungen für geeignete Prozeduren, und erklärt, welche Effekte zum Beispiel durch Strahlungen höherer Ordnung auftreten können.

#### Sehen Sie dazu auch

Wichtige Einstellungen f
ür den Betrieb (
Seite 106)

#### 6.9.1 Allgemeine Hinweise

Beachten Sie die folgenden Hinweise zur Optimierung der Geräteleistung:

- Stellen Sie sicher, dass die Qualität des Lösungsmittels mit der Anwendung kompatibel ist. In vielen Fällen sollte die Fluoreszenz-Qualität der Lösungsmittel ausreichend sein, um eine geringe Hintergrundfluoreszenz zu gewährleisten. Jedoch können ebenso normale Lösungsmittel mit HPLC-Qualität ausreichend sein.
- Im allgemeinen benötigt die Fluoreszenz-Detektion andere Lösungsmittelqualitäten im Vergleich zur UV-Detektion.
   Fluoreszierende Verunreinigungen oder Schmutzpartikel können zu Hintergrundfluoreszenz und Streulicht führen. Damit wird das Basislinienrauschen erhöht und der dynamische Bereich des Detektors verringert.
- Die Erfahrung zeigt, dass die Eignung eines Lösungsmittels für eine Anwendung stark von der gewählten Detektionswellenlänge und der benötigten Nachweisgrenze abhängt. Weiterführende Informationen und Anleitung, wie die Hintergrundfluoreszenz des Lösungsmittels bestimmt werden kann, finden Sie im Fluorescence Method Development Handbook.
- Entgasen Sie das Lösungsmittel.
- Berücksichtigen Sie die Auswirkungen der Einstellungen auf die Detektion.
Lassen Sie die Gerätetüren während des Betriebs geschlossen, um die Messzelle keinen Änderungen der Umgebungstemperatur auszusetzen, die zu einer Basisliniendrift führen können.

Behalten Sie die Betriebsdauer der Lampe im Auge, und legen Sie entsprechende Wartungsintervalle fest (siehe Überwachen der Lampenbetriebsdauer (> Seite 136)).

• Beachten Sie die allgemeinen Richtlinien und Empfehlungen zur Verwendung von Lösungsmitteln und Additiven im Chromatographie-System. Informieren Sie sich dazu in Verwenden von Lösungsmitteln und Additiven in der Vanquish-System-Betriebsanleitung.

### Sehen Sie dazu auch

- Übersicht der Parameter zur Optimierung () Seite 109)
- Überwachen der Lampenbetriebsdauer (> Seite 136)

## 6.9.2 Übersicht der Parameter zur Optimierung

Die folgende Tabelle verschafft Ihnen einen Überblick über die Parameter, die die Leistung des Detektors beeinflussen und zeigt, welche Leistungsparameter betroffen sind. Die folgenden Abschnitte enthalten Hinweise zur Auswahl der Parameter.

Parameter	Beeinflusst
Anregungswellenlänge, Emissionswellenlänge	Fluoreszenzintensität, Nachweisgrenze, Selektivität
Sensitivity	Basislinienrauschen, max. Fluoreszenzintensität
Filterrad-Einstellung	Basislinienrauschen, erlaubter Emissionswellenlängenbereich
PMT	Emissionswellenlängenbereich (und damit den Anregungswellenlängenbereich)
Ansprechzeit	Basislinienrauschen, Peakbreite, Peakhöhe
Datenaufnahmerate	Auflösung der Peaks, Speicherplatz, ggf. Basislinienrauschen
Messzellen-Temperatur	Fluoreszenzintensität, Reproduzierbarkeit
Lampen-Modus	Lebensdauer der Lampe, Basislinienrauschen, maximale Datenaufnahmerate
Basislinienverhalten	Verlauf der Basislinie nach Schalten der Wellenlänge, der Empfindlichkeit, der Filterrad- Einstellung, oder des PMT.

## 6.9.3 Optimale Anregungs- und Emissionswellenlänge bestimmen

Die wichtigsten zu optimierenden Parameter sind die Anregungs- und Emissionswellenlängen. Die folgenden Kriterien sind bei der Festlegung einer Wellenlänge für die Analyse grundsätzlich von Bedeutung:

- Wählen Sie, wenn möglich, eine Anregungswellenlänge auf dem Absorptionsmaximum der Probenbestandteile.
- Vermeiden Sie den Wellenlängenbereich, in dem Lösungsmittel absorbieren (zum Beispiel, unterhalb von 220 nm für Methanol und unterhalb von 210 nm für Acetonitril). Die Anregungswellenlänge sollte daher immer über dem UV-Cutoff des Lösungsmittels ausgewählt werden. Informationen zu den UV-Cutoff-Wellenlängen von Lösungsmitteln finden Sie unter UV-Cutoff-Wellenlängen von Lösungsmitteln (▶ Seite 190).
- Die Emissionswellenlänge muss mindestens 20 nm größer als die Anregungswellenlänge gewählt werden.

**TIPP** In Chromeleon 7 steht Ihnen das Aufzeichnen eines FL-Feldes (FL Field Acquisiton) zur Verfügung, welches die Bestimmung der Retentionszeiten und Absorptionsmaxima deutlich vereinfacht.

Weitere Informationen zur Wahl der geeigneten Wellenlängen finden Sie im mitgelieferten Fluorescence Method Development Handbook.

### Wellenlängenschaltung in der Software Chromeleon

Thermo Fisher Scientific empfiehlt, auf einem Kanal zu messen und die Wellenlängen zeitgesteuert für die nacheinander detektierten Peaks der unterschiedlichen Probenbestandteile einzustellen, anstatt im Multi-Channel-Modus alle interessanten Wellenlängen simultan zu messen.

Verwenden Sie den Instrument Method Wizard, um die Startwellenlänge für einen Kanal festzulegen. Öffnen Sie nach Beendigung des Assistenten die Methode im Instrument Method Editor. In der Modulansicht (Module View) für den Detektor können Sie auf der Registerkarte **Timetable** definieren, zu welchen Zeiten die Wellenlängen und andere Parameter geschaltet werden sollen.

Einzelheiten hierzu finden Sie auch in der Chromeleon-Hilfe.

## 6.9.4 Sensitivity (Detektorempfindlichkeit)

Mit der Empfindlichkeit kann das Signal-Rausch-Verhältnis in einem Chromatogramm optimiert werden. Je nach Fluoreszenzintensität kann es notwendig sein, die Empfindlichkeit während einer Analyse mehrfach anzupassen. Für die Wahl der optimalen Empfindlichkeit müssen die maximal auftretenden Emissionsintensitäten in einem separaten Probenlauf ermittelt werden.

## Empfindlichkeitseinstellung und Autoranging

Die Intensität des Emissionslichts nach dem Emissions-Monochromator wird mit einem oder zwei PMTs gemessen. Die Empfindlichkeit des PMT kann über die Einstellung **Sensitivity** in 8 Stufen (1 bis 8) angepasst werden. Mit jeder Stufe der Sensitivity wird die Empfindlichkeit der PMTs in etwa um den Faktor 2 größer. Hat ein Peak mit Sensitivity = 5 zum Beispiel eine Höhe von 15 Millionen Counts, dann vergrößert sich die Peakhöhe auf ca. 30 Millionen Counts, wenn mit Sensitivity = 6 gemessen wird.

- Wird die Empfindlichkeit zu klein gewählt, kann aufgrund der zu geringen Peakhöhe kein optimales Signal-Rausch-Verhältnis mehr erreicht werden (siehe Abbildung, Sensitivity = 1).
- Wird die Empfindlichkeit zu groß gewählt, dann geht das Signal des PMT in die Sättigung. Der Detektor schaltet in diesem Fall die Sensitivity automatisch zurück. Im Chromatogramm erscheinen zu Beginn und Ende des übersteuerten Bereichs Linien, die das Umschalten markieren (siehe Abbildung, Sensitivity = 8). Ist der Detektor nicht mehr übersteuert, wird die Sensitivity-Einstellung vor der Sättigung wiederhergestellt.

**TIPP** Im Chromeleon Audit Trail erscheint bei einer automatischen Reduzierung der Empfindlichkeit sowie bei der Wiederherstellung des Ursprungswertes eine Warnung, die den Wert der Empfindlichkeit nach der automatischen Anpassung angibt.



Abbildung 33: Auswirkung der Sensitivity

Beim Vanquish Fluoreszenz-Detektor wird die vom PMT gemessene Intensität des Emissionslichts auf die (von dem Referenzsensor gemessene) Bestrahlungsintensität der Probe normiert. Deshalb können die vom Emissionskanal angezeigten Werte (Counts) nicht direkt für die Optimierung der Empfindlichkeit verwendet werden. In der Abbildung erreicht der PMT bei 63 Millionen (normierten) Counts die Sättigung. Unter anderen Messbedingungen kann die Sättigung aber auch erst bei einem Wert von z.B. 100 Millionen Counts eintreten.

Verwenden Sie daher zur Optimierung der Empfindlichkeit den Parameter **MaxPMTSaturation** für jeden Peak wie unten beschrieben. Dieser zeichnet kontinuierlich die größte am PMT aufgetretene Aussteuerung in Prozent vom Sättigungswert auf seit dem letzten Chromeleon-Befehl **ClearMaxPMTSaturation**. Das Ergebnis wird aufgezeichnet in Prozent der maximal erlaubten PMT-Sättigung.

### Bestimmung der optimalen Empfindlichkeit

Bestimmen Sie die optimale Empfindlichkeit in einem separaten Probenlauf, nachdem Sie die optimalen Wellenlängen wie unten beschrieben bestimmt haben.

• Verwenden Sie dazu eine Probe (Standard) mit der maximal zu erwartenden Konzentration der Analyten.

- Wählen Sie eine Empfindlichkeit, bei welcher noch keine Übersteuerung erwartet wird (zum Beispiel 1 oder 2).
- Haben die einzelnen Peaks in einem Chromatogramm sehr unterschiedliche Höhen, dann kann das Signal-Rausch-Verhältnis der kleinen Peaks durch eine zeitgesteuerte Umschaltung der Empfindlichkeit während des Laufs zwischen den Peaks verbessert werden mit Hilfe einer Timetable.
- Nehmen Sie immer die gleiche PMT-Einstellung, um die Empfindlichkeit f
  ür die Analyse zu bestimmen.

Ein Beispiel für eine Instrumentenmethode finden Sie in der *Chromeleon-Hilfe*.

### Gehen Sie wie folgt vor

- 1. Öffnen Sie die zu optimierende Methode in dem **Script Editor**. Die notwendigen Befehle müssen manuell in der Methode eingefügt werden.
- 2. Fügen Sie einen Chromeleon-Befehl **ClearMaxPMTSaturation** zu Beginn des Laufs hinzu.
- Fügen Sie nach dem Peakmaximum eines erwarteten Peaks (oder einer Gruppe von Peaks) einen Log-Befehl für das Property MaxPMTSaturation ein, um den maximalen Sättigungswert zu diesem Zeitpunkt in den Audit Trail zu schreiben (siehe Abbildung unten).
- Fügen Sie einen ClearMaxPMTSaturation-Befehl kurz danach ein, vor der Elution des nächsten Peaks, um den MaxPMTSaturation-Wert auf null zurück zu stellen (siehe Abbildung unten).



Abbildung 34: Überwachen des Parameters MaxPMTSaturation

5. Wiederholen Sie die Schritte 3 und 4 für alle Peaks.

 Messen Sie eine Probe und lesen Sie die Werte f
ür MaxPMTSaturation in der Software Chromeleon ab und werten Sie diese aus:

Wert	Was ist zu tun
< 30%	Die Empfindlichkeit sollte erhöht werden. Als grobe Richtlinie gilt: < 30%: um eine Stufe < 15%: um zwei Stufen Wiederholen Sie dann die Optimierung ab Schritt 1.
30% - 80%	Die Empfindlichkeit ist optimal eingestellt.
80% - 99%	Die Empfindlichkeit sollte um eine Stufe reduziert werden, um eine unbeabsichtigte Übersteuerung bei Schwankungen der Konzentration zu vermeiden.
≥ 100%	Die Empfindlichkeit sollte um mindestens eine Stufe reduziert werden. In den meisten Fällen hat der Detektor die Sensitivity bereits automatisch reduziert ("Autoranging"). Sehen sie im Audit-Trail nach und wiederholen Sie mit der dort angegebenen kleinsten Empfindlichkeit nach dem letzten Autoranging die Optimierung ab Schritt 1.

Sind die maximal zu erwartenden Konzentrationen der Analyten nicht bekannt, so sollte die Empfindlichkeit bereits bei kleineren Werten von **MaxPMTSaturation** reduziert werden.

#### TIPP

- Die mit **MaxPMTSaturation** angezeigte maximale Aussteuerung bezieht sich immer auf die Signale beider PMTs bei den gegebenen Parametern aller Kanäle.
- Sie können die Fluoreszenz-Detektion komplett "ausschalten", indem Sie die Empfindlichkeit auf Off setzen. Dies kann zum Beispiel nötig sein, wenn Sie das Signal eines vorgeschalteten UV-Detektors für eine höher konzentrierte Probe aufnehmen möchten, oder einen extrem hohen Peak ausblenden möchten, der uninteressant ist aber eine automatische Anpassung der Empfindlichkeit auslöst.

## 6.9.5 Filterrad (Nur VF-Detektoren)

Damit Streulicht nicht in den PMT gelangt, werden im Lichtweg zwischen Messzelle und Emissions-Monochromator optische Kantenfilter auf einem Filterrad eingesetzt. Dabei wird Licht herausgefiltert, das durch die direkte Streuung der Lichtquelle entsteht. Zusätzlich lässt ein Gitter-Monochromator aber auch Fraktionen (die Hälfte, ein Drittel, ...) der ausgewählten Wellenlänge passieren. Wird der Emissions-Monochromator zum Beispiel auf 500 nm eingestellt, könnte auch Licht zweiter Ordnung mit 250 nm von der Probe auf den PMT gelangen. Dieses Streulicht höherer Ordnungen wird ebenfalls durch die Filter eliminiert.

Idealerweise lassen diese Filter alle Wellenlängen passieren, die größer als ihre Langpasswellenlänge sind. Licht, dessen Wellenlänge kleiner als die Langpasswellenlänge ist, wird blockiert. In der Praxis gibt es zwischen Sperr- und Durchlassbereich einen Übergangsbereich, in welchem das Licht mit zunehmender Wellenlänge immer besser transmittiert wird. Bei der Langpasswellenlänge lassen die Filter ca. 50% des Lichts passieren. Die Abbildung zeigt beispielhaft das Transmissionsverhalten des 280 nm-Filters.



Abbildung 35: Transmissionsverhalten des 280 nm-Filters

Die Filterrad-Einstellung kann automatisch oder manuell gewählt werden:

 Bei der Einstellung Auto wählt der Detektor automatisch eine Stellung des Filterrads aus. Es wird der erste Filter eingestellt, dessen Langpasswellenlänge kleiner als die eingestellte Emissionswellenlänge ist. Dieser Modus ist standardmäßig aktiviert und liefert für die meisten Anwendungen die besten Ergebnisse.

- Sie können die Filterradeinstellung (d.h., die Langpasswellenlänge des Filters) manuell wählen. Es stehen Filter mit 280 nm, 370 nm, 435 nm oder 530 nm zur Verfügung. In Sonderfällen (z.B. wenn die Emissionswellenlänge nahe an der Langpasswellenlänge eines Filters liegt) kann über die manuelle Einstellung ein besseres Ergebnis als mit der Einstellung Auto erzielt werden. Beachten Sie Folgendes:
  - Die eingestellte Emissionswellenlänge darf dabei nicht mehr als 15 nm unter der Wellenlänge des Filters liegen.
  - Bei der Einstellung Open bleibt das Filterrad geöffnet. Verwenden Sie diese Einstellung, wenn Sie mit einer Emissionswellenlänge zwischen 220 nm und 280 nm messen möchten.
  - Wählen Sie die Position des Filterrads beispielsweise in Kombination mit dem Zero-Order-Modus manuell. Hierbei wird der Emissions-Monochromator geöffnet und das emittierte Licht über den gesamten Wellenlängenbereich gemessen. Nutzen Sie den Langpassfilter dazu, Licht unterhalb des Wellenlängenbereichs auszublenden, der für die Anwendung interessant ist. Beachten Sie bei dieser Einstellung, dass diese Einstellung üblicherweise aufgrund des Streulichts und der Autofluoreszenz des Eluenten eine geringere Empfindlichkeit benötigt (der gesamte spektrale Bereich wird aufgezeichnet).

Weitere Informationen zur Filterrad-Einstellung finden Sie im mitgelieferten Fluorescence Method Development Handbook.

## 6.9.6 PMT (nur wenn zweiter PMT installiert ist)

Die Detektoren können mit einem zweiten PMT für den nahen Infrarotbereich (bis 900 nm) ausgestattet sein. Wird die PMT-Einstellung auf **Auto** gesetzt, wählt der Detektor den geeigneten PMT für jede Messung aus. Alternativ können Sie manuell wählen, welcher PMT verwendet werden sollte. Verwenden Sie PMT1 für Messungen im UV- / VIS-Bereich. Verwenden Sie PMT2 für Messungen im Infrarot-Bereich.

**TIPP** Kalibrierung und Quantifizierung müssen immer mit derselben PMT-Einstellung durchgeführt werden. Eine Kalibrierung mit PMT1 ist für Messungen mit PMT2 nicht zwingend gültig.

## 6.9.7 Datenaufnahmerate und Ansprechzeit (Single-Channel-Modus)

Die Software Chromeleon berechnet automatisch die beste Ansprechzeit (Response Time), basierend auf dem Wert, den Sie für die Datenaufnahmerate im Instrument Method Wizard eingeben. Beachten Sie die folgenden Hinweise zur Auswahl der Datenaufnahmerate. Wenn Sie eine abweichende Ansprechzeit wählen möchten, finden Sie im folgenden Abschnitt ebenfalls Hinweise dazu.

**TIPP** Im Multi-Channel-Modus werden die Ansprechzeit und Datenaufnahmerate immer automatisch festgelegt. Sie können diese Parameter beeinflussen, indem Sie eine andere Messqualität (Multi-Channel Performance) wählen.

## Datenaufnahmerate

Die Datenaufnahmerate gibt an, wie viele Datenpunkte pro Sekunde (Hz) die Software Chromeleon vom Detektor übernimmt und als Rohdaten abspeichert.

Wenn Sie eine Datenaufnahmerate wählen, beachten Sie folgende Hinweise:

- Jeder Peak sollte in der Regel durch wenigstens 20 Datenpunkte definiert werden. Für Chromatogramme mit koeluierenden Peaks oder einem geringen Signal/Rausch-Verhältnis werden 40 Datenpunkte pro Peak empfohlen.
- Bei einer zu niedrigen Datenaufnahmerate werden Peakanfang, Maxima und Peakende nicht genau bestimmt.
- Wenn die Datenaufnahmerate zu hoch ist, benötigen die Daten gegebenenfalls mehr Speicherplatz als nötig und die Verarbeitungszeit für die Post-Run-Analyse verlängert sich.
- Wählen Sie eine niedrigere Datenaufnahmerate (beispielsweise 1,0 Hz), wenn alle Peaks verhältnismäßig breit sind. Dies spart Speicherplatz und beschleunigt die Darstellung der Daten in der Software Chromeleon.
- Wählen Sie eine höhere Datenaufnahmerate (beispielsweise 10,0 Hz), wenn die interessanten Peaks weniger als einige Sekunden breit sind.
- Der gewählte Lampen-Modus hat ebenfalls Auswirkungen auf die maximale Datenaufnahmerate.

### Ansprechzeit

Die Ansprechzeit ist ein Maß dafür, wie schnell der Detektor auf eine Signaländerung reagiert.

- Wählen Sie eine Ansprechzeit, die bei circa 1/3 der Peakbreite bei halber Peakhöhe des schmalsten interessanten Peaks liegt. Bei einer längeren Ansprechzeit können mehr Signale gemittelt werden, wodurch das Kurzzeitrauschen reduziert und das Signal-Rausch-Verhältnis verbessert wird.
- Wenn eine zu lange Ansprechzeit gewählt wird, kann dies zu reduzierten Peakhöhen und asymmetrischen Peakformen führen. Dicht hintereinander eluierende Peaks können ggf. nicht mehr korrekt getrennt werden. Eine korrekt gewählte Ansprechzeit führt zu einer wesentlichen Reduzierung des Basislinienrauschens, reduziert die Peakhöhe jedoch nur geringfügig.

**TIPP** Die beste Kombination von Datenaufnahmerate, Ansprechzeit und Peakbreite erhalten Sie, wenn Sie das Kontrollkästchen **Link data collection parameters** in den Detektoreigenschaften des Chromeleon Instrument Method Wizards oder Chromeleon Instrument Method Editors auswählen.

#### Sehen Sie dazu auch

■ Lampen-Modus (▶ Seite 119)

## 6.9.8 Messzellen-Temperatur

Die Temperatur der Probe hat Einfluss auf deren Fluoreszenz. Bei Fluoreszenz-Detektion besteht daher das Problem, dass Messergebnisse durch Schwankungen der Umgebungstemperatur stark beeinflusst werden können. Vanquish Fluoreszenz-Detektoren bieten die Möglichkeit, die Messzelle zu heizen und somit die Temperatur in der Messzelle konstant zu halten, selbst wenn die Raumtemperatur variiert.

Die Fluoreszenz der meisten Substanzen nimmt mit steigender Temperatur ab. Deshalb sollte die Temperatur nicht zu hoch ausgewählt werden. Sie muss jedoch über der Temperatur liegen, die in der Optik des Detektors vorherrscht, die wiederum von der Umgebungstemperatur beeinflusst wird. Wählen Sie die Temperatur daher etwa 15°C über der zu erwartenden Umgebungstemperatur.

*Beispiel*: Im Labor wird eine maximale Temperatur von 27°C erwartet. Setzen Sie die Messzellen-Temperatur auf 42°C.

**TIPP** Thermo Fisher Scientific empfiehlt, die Messzellen-Temperatur über den Signalkanal **FLD\_FlowCell** immer mit aufzuzeichnen. Im Fall einer Störung kann das Temperatursignal hilfreiche Hinweise auf die Ursache der Störung liefern.

## 6.9.9 Lampen-Modus

Für die Xenon-Blitzlampe des Detektors stehen drei unterschiedliche Modi zur Verfügung. Durch die Wahl eines anderen Lampenmodus in Phasen, in denen keine interessanten Peaks eluieren, kann die Lebensdauer der Lampe verlängert werden.

Bei der Berechnung der Lebensdauer der Lampe wird nur die Zeit mit einberechnet, in der die Lampe auch blitzt. Dies entspricht ungefähr der Laufzeit der Datenaufnahme.

## Verlängerung der Lebensdauer der Lampe

Sie können Folgendes tun, um die Lebensdauer der Lampe zu verlängern:

- Schalten Sie die Lampe aus, indem Sie die Datenaufnahme beenden, wenn die Basislinie für Sie nicht mehr interessant ist. Es ist im Allgemeinen nicht erforderlich, die Lampe während der gesamten chromatographischen Trennung eingeschaltet zu lassen.
- Schalten Sie während der chromatographischen Trennung zwischen verschiedenen Lampen-Modi um, zum Beispiel vom Modus LongLife (zwischen den Peaks) in den Modus Standard oder HighPower (für die interessanten Peaks). Dadurch kann die Lebensdauer der Lampe erhöht werden, ohne dass die Messung an den wichtigen Stellen an Empfindlichkeit einbüßt.

Verwenden Sie den Instrument Method Wizard, um den Basis-Lampenmodus festzulegen. Öffnen Sie nach Beendigung des Assistenten die Methode im Instrument Method Editor. In der Modulansicht (Module View) für den Detektor können Sie auf der Registerkarte **Timetable** definieren, zu welchen Zeiten der Lampen-Modus sowie andere Parameter geschaltet werden sollen.

Ein Beispiel zur geeigneten Wahl des Lampen-Modus finden Sie im Fluorescence Method Development Handbook.

### HighPower-Modus

Die Lampe blitzt mit der höchstmöglichen Frequenz von 300 Hz. Die Empfindlichkeit ist etwa doppelt so hoch wie im Modus **Standard**. Die zu erwartende Lebensdauer der Lampe beträgt etwa 1300 Stunden.

Dieser Modus wird für Anwendungen empfohlen, für die eine höchste Empfindlichkeit gewünscht wird.

Modus Standard	
	Die Lampe blitzt mit einer mittleren Frequenz von 100 Hz. Die zu erwartende Lebensdauer der Lampe beträgt etwa 4000 Stunden. Dieser Modus unterstützt Datenaufnahmeraten bis 100 Hz.
	Dieser Modus wird für Anwendungen empfohlen, für die eine hohe Empfindlichkeit gewünscht wird.
Modus LongLife	
	Die Lampe blitzt mit einer langsameren Frequenz von 20 Hz. Wenn Sie in diesem Lampen-Modus arbeiten, können Sie die Lebensdauer einer Lampe gegenüber dem Modus <b>Standard</b> in etwa vervierfachen. Die zu erwartende Lebensdauer der Lampe beträgt etwa 16000 Stunden. Beachten Sie aber, dass sich das Signal-Rausch-Verhältnis im Vergleich zum Modus <b>Standard</b> etwa auf die Hälfte reduziert. Dieser Modus unterstützt nur Datenaufnahmeraten bis 20 Hz.
	Dieser Modus wird für Bereiche des Chromatogramms empfohlen, in denen das Signal-Rausch-Verhältnis eine geringere Rolle spielt, zum Beispiel:

- Wenn keine interessanten Peaks eluieren
- Während der Waschzyklen und Reäquilibrierzeit

## 6.9.10 Basislinienverhalten

Werden im Laufe einer Methode Wellenlängen geschaltet oder auch die Empfindlichkeit, Filterrad-Einstellung oder die PMT-Einstellung verändert, so kann es im Chromatogramm aufgrund der Hintergrundfluoreszenz bei der neuen Wellenlängeneinstellung zu Sprüngen im Basislinienverlauf kommen. Drei verschiedene Basislinien-Modus-Einstellungen definieren wie die Basislinie sich in diesen Fällen verhält:

- Zero: Die Basislinie wird auf Null gesetzt
- Append: Die Basislinie wird an das bisherige Signal angehängt
- Free: Die Basislinie "springt" auf den aktuellen absoluten Wert

**TIPP** Sie können jederzeit über den Befehl **AutoZero** einen Nullabgleich durchführen, d.h. die gemessene Intensität zum Zeitpunkt des Autozero wird von den in der Folge gemessenen Intensitäten subtrahiert. Wählen Sie für den Nullabgleich immer einen Zeitpunkt, in dem keine Probe durch die Messzelle fließt. Die Werte werden gespeichert und können über den Befehl **ClearAutoZero** wieder zurückgesetzt werden.

## 6.10 Außerbetriebnahme des Gerätes

Wird das Gerät einige Zeit lang nicht betrieben, folgen Sie den Anweisungen in diesem Abschnitt, um das Gerät außer Betrieb zu nehmen.

**TIPP** Mit der Software Chromeleon können bestimmte Abläufe zur Betriebsunterbrechung des Chromatographie-Systems automatisch durchgeführt werden. Dies gilt zum Beispiel für das Herunterfahren der Flussrate, Herabsetzen der Temperatur in temperaturgesteuerten Geräten und das Ausschalten der Detektorlampen. Weitere Informationen zum **Smart Shutdown** und **Smart Standby** finden Sie in der *Chromeleon-Hilfe*.

## 6.10.1 Kurzzeitige Außerbetriebnahme (Betriebsunterbrechung)

Um den Gerätebetrieb für kurze Zeit zu unterbrechen (kurzzeitige Außerbetriebnahme), zum Beispiel über Nacht, beachten Sie, abhängig vom Systemaufbau, die folgenden Hinweise für die Vanquish-Systemmodule:

Detektortyp	Beschreibung
Charged-Aerosol- Detektor	Stellen Sie sicher, dass ausreichend Gas zur Verfügung steht, um den Gasfluss durch den Detektor eingeschaltet zu lassen. Dies verhindert, dass sich Lösungsmittel- oder Probenrückstände im Detektor ansammeln. Der Gasfluss muss eingeschaltet sein, wenn der Pumpenfluss zum Detektor eingeschaltet ist.
UV/VIS-Detektoren:	Die Lampe (bzw. Lampen) im Detektor können eingeschaltet bleiben. Nur Variabler Wellenlängendetektor und VH-D10 Diodenarray-Detektor: Der Shutter kann zum Schutz der Messzelle geschlossen werden.
Fluoreszenz- Detektor	Schalten Sie die Temperatursteuerung für die Messzelle aus.

• Beachten Sie Folgendes für Ihren Vanquish-Detektor:

• Fördern Sie ein geeignetes Lösungsmittel mit einer Flussrate von 0,05 mL/min.

Prüfen Sie die untere Druckgrenze für die Pumpe und passen Sie den Wert gegebenenfalls an. Wenn der Druck auf einen Wert unterhalb des unteren Grenzwertes fällt, schaltet die Pumpe den Fluss ab.

• Stellen Sie das Injektionsventil im Autosampler auf die Position Inject.

- Stellen Sie sicher, dass die S\u00e4ulentemperatur nicht mehr als 40 °C betr\u00e4gt.
- Warten Sie bei Wiederaufnahme des Betriebs, bis sich der Fluss äquilibriert hat, und vergewissern Sie sich, dass die Betriebsparameter der anderen Module auf geeignete Werte eingestellt sind, ehe Sie fortfahren.

## 6.10.2 Langfristige Außerbetriebnahme

#### Außerbetriebnahme des Gerätes

Folgen Sie bei längeren Betriebsunterbrechungen den folgenden Anweisungen.

**TIPP** Die Außerbetriebnahme des Geräts beeinflusst den Betrieb Ihres Systems. Wenn Sie das Gerät außer Betrieb nehmen, beachten Sie auch die Anweisungen für die Außerbetriebnahme anderer Vanquish-Systemmodule und führen Sie die entsprechenden Schritte durch (siehe dazu die *Betriebsanleitungen* für die Module).

- 1. Schalten Sie den Pumpenfluss ab.
- 2. Entfernen Sie die Säule.
- Verbinden Sie die freien Enden der Kapillaren des Säulenthermostaten mit einem Verbindungsstück (zum Beispiel, dem Viper-Verbindungsstück aus dem Zubehörkit des Systems) und starten Sie den Pumpenfluss wieder.
- 4. Spülen Sie die Messzelle mit einem geeigneten Lösungsmittel (mindestens HPLC-Qualität). Beachten Sie Folgendes:

**TIPP** Bei einem Vanquish Core-System, das zur Verwendung für Normal-Phasen-kompatible Lösungsmittel und Additive modifiziert ist, beachten Sie die Informationen über die Spülflüssigkeit im Kapitel *Besonderheiten bei Normal-Phase kompatiblen Lösungsmitteln und Additiven* in der *Betriebsanleitung* für die Pumpe.

Situation nach der Außerbetriebnahme	Wenn kein Additiv verwendet wird	Wenn ein Additiv verwendet wird
Gerät und Messzelle bleiben nach der Außerbetriebnahme im Labor	Spülen Sie das System zum Beispiel mit Methanol. 100%-iges Acetonitril sollte nicht verwendet werden.	Spülen Sie das System einige Male (zum Beispiel mit 1,0 mL/min 10 Minuten lang bei einem Standard-System) mit einem Gemisch aus Methanol und Wasser (50:50), um eine Aufkonzentrierung von Salzen im Flusspfad zu vermeiden. Wenn die Lösungsmittel in der Messzelle nicht mit Wasser mischbar sind, verwenden Sie einen geeigneten Löslichkeitsvermittler.
Das Gerät und die Messzelle sollen nach der Außerbetriebnahme transportiert oder verschickt werden	Spülen Sie das System mit Isopropanol.	Spülen Sie das System zunächst einige Male (zum Beispiel mit 1,0 mL/ min 10 Minuten lang bei einem Standard-System) mit einem Gemisch aus Methanol und Wasser (50:50), um eine Aufkonzentrierung von Salzen im Flusspfad zu vermeiden. Wenn die Lösungsmittel in der Messzelle nicht mit Wasser mischbar sind, verwenden Sie einen geeigneten Löslichkeitsvermittler. Spülen Sie anschließend das System mit Isopropanol.

#### ACHTUNG

Probenreste, Verunreinigungen von der Säule oder Puffer mit hohen Salzkonzentrationen können sich in der Messzelle ablagern. Dies kann zu Schäden an der Messzelle führen. Außerdem können säurehaltige Lösungsmittel die Messzelle beschädigen.

- Spülen Sie die Messzelle mit einem geeigneten Lösungsmittel, bevor Sie den Betrieb unterbrechen.
- Füllen Sie die Messzelle mit Hilfe des Spül- und Injektionskits mit Isopropanol.
- Die Messzelle sollte nicht mit reinem Wasser gefüllt sein, um Algenbildung zu vermeiden. Wenn Sie die Messzelle mit Wasser füllen möchten, mischen Sie dem Wasser 10% Isopropanol in HPLC-Qualität bei.
- 5. Schalten Sie den Pumpenfluss zum Gerät ab. Warten Sie, bis das System druckfrei ist, bevor Sie mit der Außerbetriebnahme des Geräts fortfahren.
- 6. Entfernen Sie die Kapillaren vom Einlass und Auslass der Messzelle.
- Verschließen Sie nach dem Entfernen der Flussverbindungen den Einlass und den Auslass der Messzelle mit den Verschlussstopfen, mit denen die Messzelle ausgeliefert wurde. Schützen Sie die Einlasskapillaren mit Verschlusskappen.
- 8. Dieser Schritt ist wie folgt abhängig:

Situation	Schritte
Das Gerät und alle anderen Systemmodule bleiben im Systemturm und alle Systemmodule sollen ausgeschaltet werden	Schalten Sie das System über den System-Ein-/Ausschalter am Systemsockel aus.
Das Gerät soll nach der Außerbetriebnahme transportiert oder verschickt werden	Wenn eines der Module aus dem Systemturm entfernt werden soll, schalten Sie <i>alle</i> Systemmodule über ihren Hauptschalter aus. Das Drücken des System-Ein- /Ausschalters reicht nicht aus, um die Geräte vollständig auszuschalten.
	Folgen Sie den Anweisungen in Transportieren und Versenden des Gerätes (▶ Seite 152).

**TIPP** Um bei Betriebsunterbrechungen Schäden an der Detektoroptik durch Staubpartikel zu vermeiden, achten Sie darauf, dass am Detektor immer eine Messzelle oder die mitgelieferte Messzellen-Abdeckung installiert ist.

#### Sehen Sie dazu auch

- Anschließen der Einlasskapillare (> Seite 72)
- Anschließen der Detektor-Waste-Leitung (> Seite 76)

## 6.10.3 Wiederaufnehmen des Gerätebetriebs nach langfristiger Außerbetriebnahme

Um das Gerät nach einer langfristigen Außerbetriebnahme wieder in Betrieb zu nehmen, gehen Sie wie folgt vor:

- 1. Bereiten Sie die anderen Module im Vanquish-System vor und starten Sie diese; folgen Sie dabei den Anweisungen in den *Betriebsanleitungen* für die Module. Beachten Sie besonders den Abschnitt *Vorbereiten des Moduls für den Betrieb*.
- 2. Spülen Sie die Komponenten, die im Flussweg vor der Messzelle angeschlossen sind, bevor Sie die Flussverbindungen des Systems zur Messzelle herstellen.
- 3. Verbinden Sie die Einlasskapillare mit dem Einlass der Messzelle und die Waste-Leitung mit dem Auslass der Messzelle.
- 4. Schalten Sie das Gerät ein. Beachten Sie Folgendes:

Situation	Aktion
Wenn das Gerät im Systemturm	Schalten Sie das System über den
geblieben ist und alle Systemmodule	System-Ein-/Ausschalter am
abgeschaltet wurden.	Systemsockel ein.
Wenn der Gerätebetrieb nach einem	Schalten Sie das Gerät über den
Transport wieder aufgenommen wird.	Hauptnetzschalter ein.

5. Bevor Sie mit einer Analyse beginnen, warten Sie, bis sich der Detektor äquilibriert hat, und vergewissern Sie sich, dass er für den Betrieb bereit ist.

# 7 Wartung und Service

Dieses Kapitel enthält Informationen zur routinemäßigen Wartung und zu Servicearbeiten, die Sie als Anwender durchführen können.

## 7.1 Einführung in Wartung und Service

Dieses Kapitel enthält Informationen zur routinemäßigen Wartung sowie zu Service- und Reparaturarbeiten, die Sie als Anwender durchführen können.



Weiterführende Wartungs- oder Servicearbeiten dürfen nur von Service-Personal durchgeführt werden, das von Thermo Fisher Scientific entsprechend zertifiziert wurde (im Folgenden kurz als Thermo Fisher Scientific-Servicetechniker bezeichnet).

Das Gerät ist für einfache Wartungs- und Reparaturarbeiten ausgelegt. Die Komponenten, für die der Anwender Servicearbeiten durchführen kann, sind von der Vorderseite aus zugänglich. Wenn nicht anders angegeben, kann das Gerät für die Wartungsarbeiten im System bleiben.

Es ist nicht erforderlich, die Türen für die Wartungsarbeiten zu entfernen. Sie können jedoch eine Tür entfernen, sollte dies aus einem bestimmten Grund oder für bestimmte Arbeiten erforderlich sein. Wenn Sie eine Tür entfernen müssen, folgen Sie den entsprechenden Schritten in Tauschen der Türen (▶ Seite 150).

## 7.2 Sicherheitshinweise zu Wartung und Service

Beachten Sie bei der Durchführung von Wartungs- und Servicearbeiten folgende Sicherheitshinweise:



Beachten Sie alle Warnhinweise und Vorsichtsmaßnahmen in Sicherheitsmaßnahmen (▶ Seite 21).



## WARNUNG—Hochspannung

Im Gerät treten hohe Spannungen auf, die zu einem Stromschlag führen können.

Öffnen Sie nicht das Gehäuse oder entfernen Sie keine Schutzabdeckungen, es sei denn, Sie werden in dieser Anleitung dazu ausdrücklich aufgefordert.



## WARNUNG—Austreten gefährlicher Substanzen aus Flussverbindungen

Fluss- und Kapillarverbindungen können mit Substanzen gefüllt sein, die ein Gesundheitsrisiko darstellen können. Lösungsmittel können herausspritzen, wenn Kapillaren bersten, aus ihren Fittingen rutschen, oder nicht korrekt festgezogen sind, oder wenn Kapillarverbindungen aus anderen Gründen offen sind.

- Tragen Sie eine geeignete Schutzausrüstung und folgen Sie der Guten Laborpraxis.
- Spülen Sie schädliche Substanzen vor Beginn der Wartungs- oder Servicearbeiten mit einem geeigneten Lösungsmittel aus.



#### WARNUNG—Kippende Flüssigkeitsbehälter

Die Behälter im Solvent Rack können Flüssigkeiten mit schädlichen Substanzen enthalten. Ein Verschütten dieser Substanzen kann Gesundheits- und Sicherheitsrisiken darstellen.

Um zu vermeiden, dass die Behälter kippen, ziehen Sie bei Wartungsarbeiten nicht an den Flüssigkeitsleitungen.



#### VORSICHT—Austretende Lösungsmittel

Lösungsmittel können austreten, wenn sie unter hohem Druck stehen.

- Schalten Sie den Pumpenfluss ab, bevor Sie den Flussweg öffnen.
- Warten Sie, bis das System druckfrei ist.
- Tragen Sie eine geeignete Schutzausrüstung, wenn Sie Verbindungen im Flussweg öffnen.



#### VORSICHT—Heiße Oberflächen

Oberflächen im Inneren Geräts können sich während des Betriebs erhitzen. Das Berühren von heißen Komponenten kann zu Verbrennungen führen.

Warten Sie, bis die heißen Oberflächen abgekühlt sind, bevor Sie mit Austausch- und Wartungsarbeiten beginnen.



#### VORSICHT—Hydrostatischer Druck

Lösungsmittel können austreten, wenn Sie Verbindungen im Flussweg öffnen. Grund dafür ist der hydrostatische Druck im System, wenn Lösungsmittelbehälter oberhalb des Pumpenausgangs stehen. Bevor Sie eine Verbindung im Flussweg öffnen:

- Schalten Sie den Pumpenfluss ab und warten Sie, bis das System druckfrei ist.
- Schrauben Sie die Deckel der Lösungsmittelbehälter ab und entfernen Sie die Lösungsmittelschläuche zusammen mit den Deckeln aus den Behältern.
- Entleeren Sie die Lösungsmittelschläuche. Einzelheiten entnehmen Sie bitte der *Betriebsanleitung* für die Pumpe.
- Drehen Sie die Deckel der Behälter wieder fest.



#### VORSICHT—Stromschlag oder Schäden am Gerät

Auch im ausgeschalteten Zustand fließt im Gerät Strom, solange das Netzkabel eingesteckt ist. Reparaturen am Gerät, während das Gerät am Stromnetz angeschlossen ist, können zu Personenschäden führen.

- Ziehen Sie immer das Netzkabel ab, bevor Sie Reparaturen im Geräteinneren durchführen.
- Sollten Sie Gehäuseabdeckungen oder Seitenwände entfernen müssen, schließen Sie das Netzkabel keinesfalls an das Gerät an, solange die Abdeckungen und Seitenwände noch nicht montiert sind.

## 7.3 Allgemeine Regeln für Wartung und Service

Damit die Wartungs- und Servicearbeiten erfolgreich sind, beachten Sie folgende Regeln und Empfehlungen:

- Bevor Sie Service- oder Wartungsarbeiten ausführen, nehmen Sie das Gerät außer Betrieb, wenn Sie dazu aufgefordert werden.
- Verwenden Sie ausschließlich die Ersatzteile, die von Thermo Fisher Scientific ausdrücklich für das Gerät autorisiert und freigegeben sind.
- Folgen Sie allen Anweisungen Schritt für Schritt und verwenden Sie die in der Anleitung empfohlenen Werkzeuge.
- Bevor Sie eine Verbindung im Flussweg öffnen, um Kapillaren im System auszutauschen, schalten Sie den Pumpenfluss ab und warten Sie, bis das System druckfrei ist.
- Verunreinigte Komponenten können zu einer Verunreinigung des Chromatographie-Systems führen. Verunreinigungen führen zu einer schlechten Leistung der Module und des gesamten Systems oder sogar zu Schäden an den Modulen und dem System. Daher gilt:
  - Tragen Sie immer geeignete Schutzhandschuhe.
  - Legen Sie die Komponenten nur auf einer sauberen, fusselfreien Arbeitsfläche ab.
  - Halten Sie die Werkzeuge sauber.
  - Verwenden Sie zur Reinigung nur ein fusselfreies Tuch.
- Wenn Sie das Gerät zur Reparatur zurückschicken müssen, folgen Sie den Anweisungen in Transportieren und Versenden des Gerätes (> Seite 152).

## Sehen Sie dazu auch

Ersatzteile und Verbrauchsmaterialien ( Seite 185)

## 7.4 Wartung und Wartungsintervalle

Optimale Leistungsfähigkeit und maximale Verfügbarkeit des Gerätes sowie zuverlässige Ergebnisse können nur sichergestellt werden, wenn das Gerät in einem guten Zustand ist und ordnungsgemäß gewartet wird.

## 7.4.1 Wartungszeitplan

Führen Sie regelmäßig die Wartungsarbeiten in der Tabelle durch. Die in der Tabelle angegebene Häufigkeit dient als Orientierungshilfe. Welche Zeitabstände für Wartungsarbeiten optimal sind, hängt von mehreren Faktoren ab, wie den Arten und Mengen der Proben und Lösungsmittel, die mit dem Gerät verwendet werden.

Häufigkeit	Was ist zu tun
Täglich	<ul> <li>Prüfen Sie die Kapillar- und Schlauchverbindungen auf Anzeichen von Undichtigkeiten oder Blockierung.</li> </ul>
	<ul> <li>Wenn Sie mit Puffern oder Salzlösungen arbeiten, spülen Sie das Gerät nach Abschluss der Arbeiten gründlich mit einem geeigneten Lösungsmittel, das keine Puffer oder Salze enthält. Achten Sie auf den Gegendruck an der Messzelle.</li> </ul>
Regelmäßig	<ul> <li>Prüfen Sie die Flussverbindungen auf Anzeichen von Beschädigungen, wie Knicke, Risse, Schnitte oder Blockierung.</li> </ul>
	<ul> <li>Führen Sie eine Qualifizierung durch und überprüfen Sie die Betriebsdauer der Lampe je nach Erfordernis der jeweiligen Anwendung.</li> </ul>
	<ul> <li>Vergewissern Sie sich, dass auf dem Gerät noch alle Warnaufkleber vorhanden und deutlich lesbar sind.</li> <li>Sollte dies nicht der Fall sein, wenden Sie sich für Ersatzaufkleber an Thermo Fisher Scientific.</li> </ul>
Jährlich	Lassen Sie ein Mal pro Jahr vorbeugende Wartungsmaßnahmen von einem Thermo Fisher Scientific- Servicetechniker durchführen.

**TIPP** Die Software Chromeleon unterstützt Funktionen zur Überwachung und Aufzeichnung von Informationen zu Service- und Qualifizierungsarbeiten (siehe Überwachen der Lampenbetriebsdauer

( Seite 136)).

## 7.4.2 Reinigen oder Dekontaminieren des Gerätes

Das Reinigen und Dekontaminieren muss immer von qualifiziertem Personal in geeigneter Schutzausrüstung durchgeführt werden. Beachten Sie stets landesspezifische und lokale Bestimmungen.

## ACHTUNG

Wischen Sie alle auf dem System verschütteten Flüssigkeiten sofort auf. Eine längere Einwirkung kann Schäden verursachen.

## Dekontaminierung

Eine Dekontaminierung ist zum Beispiel erforderlich, wenn eine Undichtigkeit aufgetreten ist oder Flüssigkeit verschüttet wurde, oder vor Wartung oder Transport des Gerätes. Verwenden Sie ein geeignetes Reinigungs- oder Desinfektionsmittel, um sicherzustellen, dass das Gerät nach der Behandlung sicher gehandhabt werden kann.

## Erforderliche Teile

- Geeignetes Reinigungsmittel (oder Desinfektionsmittel)
- Gereinigtes Wasser
- Fusselfreie Tücher oder Papiertücher



## VORSICHT—Explosive Gasmischungen aus alkoholhaltigen Reinigungsmitteln

Alkoholhaltige Reinigungsmittel können an der Luft entzündliche und explosive Gasmischungen erzeugen.

- Verwenden Sie solche Reinigungsmittel nur, wenn erforderlich und nur in ausreichend belüfteten Räumen.
- Vermeiden Sie offene Flammen oder übermäßige Wärmeeinwirkung während der Reinigung.
- Wischen Sie die gereinigten Komponenten nach der Reinigung gründlich trocken. Betreiben Sie das Gerät nicht, wenn es nicht vollständig trocken ist.

#### ACHTUNG

Beachten Sie Folgendes:

- Verwenden Sie nur Reinigungsmittel, welche die Systemoberflächen nicht beschädigen.
- Verwenden Sie zur Reinigung der Oberflächen niemals scharfe Werkzeuge oder Bürsten.
- Verwenden Sie keine Sprays für die Reinigung.
- Achten Sie darauf, dass kein Reinigungsmittel in den Flussweg gelangt.
- Verwenden Sie kein übermäßig nasses Tuch oder Feuchttücher für die Reinigung. Achten Sie darauf, dass keine Flüssigkeiten in die funktionalen Bauteile des Geräts gelangen. Flüssigkeiten können einen Kurzschluss auslösen, wenn sie in Kontakt mit elektronischen Komponenten kommen.

#### ACHTUNG—Messzellenschacht

Die optischen Ports und der Kontaktpunkt für den Identifikationschip im Messzellenschacht sind empfindlich gegen elektrostatische Entladungen, Verunreinigungen und Kratzer.

Berühren Sie keinesfalls die optischen Ports oder andere Flächen im Messzellenschacht.

### Vorbereitungen

1. Schalten Sie das Gerät aus und trennen Sie das Netzkabel von der Stromversorgung.

#### Gehen Sie wie folgt vor

- Wischen Sie die Oberflächen mit einem sauberen, trockenen, weichen, fusselfreien Tuch oder Reinigungstuch ab. Feuchten Sie das Tuch oder Reinigungstuch gegebenenfalls mit einer Lösung aus lauwarmem Wasser und einem geeigneten Reinigungsmittel an.
- 2. Lassen Sie das Reinigungsmittel wie vom Hersteller empfohlen einwirken.
- 3. Wischen Sie alle gereinigten Oberflächen mit gereinigtem Wasser nach, damit alle Reinigungsmittelreste entfernt werden.
- 4. Trocknen Sie die Oberflächen mit einem weichen, fusselfreien Tuch oder Reinigungstuch.

## 7.4.3 Predictive Performance

Die Software Chromeleon unterstützt Funktionen zur Überwachung und Aufzeichnung von Informationen zu Service- und Qualifizierungsarbeiten am Gerät.

## Überwachen von Qualifizierungs- und Service-Intervallen

Auf speziellen Panels zu Service und Qualifizierung können Sie Intervalle für Servicearbeiten oder Qualifizierungsvorgänge festlegen. Mit diesen Funktionen, die als Predictive Performance bezeichnet werden, können Sie diese Arbeiten basierend auf den tatsächlichen Betriebs- und Nutzungsbedingungen des Gerätes planen. Sie können darüber hinaus Grenzwerte für Warnungen festlegen, um vor oder bei Ablauf eines Termins für eine Servicearbeit oder Qualifizierung daran erinnert zu werden.

Auf speziellen Panels zeigen farblich gekennzeichnete Balken die Informationen zum Zustand optisch an, damit Sie den Zustand leicht prüfen und überwachen können. Wurde ein Grenzwert für eine Warnung festgelegt, macht Sie eine Meldung im Chromeleon Audit Trail auf die Fälligkeit aufmerksam.

Service- und Qualifizierungszähler können auf null zurück gestellt werden, nachdem die erforderliche Tätigkeit ausgeführt wurde. Damit die Informationen zur Predictive Performance aktuell sind, denken Sie daran, die Zähler zurück zu stellen, nachdem Sie eine Servicearbeit oder Qualifizierung durchgeführt haben.

Weitere Informationen finden Sie in der Chromeleon-Hilfe.

## 7.4.4 Überwachen der Lampenbetriebsdauer

Die Software Chromeleon unterstützt Funktionen zur Überwachung des Lampenalters. Diese Funktion kann dabei helfen festzustellen, wann ein Lampentausch fällig ist.

Die Blitzlampe kann etwa 1-1,5x10<sup>9</sup> Blitze produzieren, bevor sie getauscht werden muss und auch nur dann, wenn die Leistung des Detektors nicht mehr ausreichend ist. Deshalb ist die Lebensdauer der Blitzlampe von Blitzfrequenz und Lampen-Modus abhängig. Anhand der Anzahl der Lampenblitze berechnet die Software Chromeleon einen Wert, der angibt, wie viele Stunden die Lampe betrieben wurde. Dabei geht die Berechnung vom Lampen-Modus **Standard** aus. Für die Berechnung der Betriebsdauer wird eine Stunde unterschiedlich gewichtet, abhängig vom gewählten Lampen-Modus.

Modus	1 Betriebsstunde wird gezählt als	Zu erwartende Lebensdauer
LongLife	0,25 Stunden	circa 16000 Stunden
Standard	1 Stunde	circa 4000 Stunden
HighPower	3 Stunden	circa 1300 Stunden

*Beispiele*: Wenn Chromeleon für die Betriebsdauer der Lampe 2000 Stunden angibt, bedeutet dies, dass die Lampe etwa die Hälfte ihrer zu erwartenden Lebensdauer erreicht hat. Wenn Sie die Lampe im Modus **LongLife** betreiben, sind also in etwa noch weitere 8000 Betriebsstunden zu erwarten.

Wählen Sie in Chromeleon den Parameter **EffectiveLampAge**. Wenn der Wert 4000 Stunden überschritten hat, sollte die Lampe ausgetauscht werden. Beachten Sie, dass der Tausch der Lampe nur von einem Thermo Fisher Scientific-Servicetechniker durchgeführt werden darf. Der Servicetechniker wird den Zähler für das Lampenalter auf null zurücksetzen, nachdem er die Lampe ausgetauscht hat.

**TIPP** Sie können die Lebensdauer der Xenon-Blitzlampe verlängern, wenn Sie während einer chromatographischen Analyse zwischen den Lampen-Modi umschalten.

## Sehen Sie dazu auch

Lampen-Modus ( Seite 119)

## 7.5 Durchführung einer Wellenlängen-Kalibrierung

Ein einfacher Initialisierungsvorgang wird nach dem Einschalten des Detektors durchgeführt. Hierfür muss eine Messzelle im Detektor installiert sein. Einzelheiten finden Sie unter Ein- und Ausschalten (▶ Seite 97).

Um eine optimale Funktionsfähigkeit und Wellenlängengenauigkeit zu gewährleisten, führen Sie die Wellenlängenkalibrierung mit Wasser (Raman-Messung) durch. Bei der Wellenlängenkalibrierung wird das Lichtspektrum der Xenon-Blitzlampe zur Kalibrierung der Anregungswellenlänge verwendet. Anschließend wird der Emissions-Monochromator anhand des Maximums des Raman-Emissionspektrums (397 nm) justiert.

## Wann

- Nach einer Installation oder einem Austausch der Messzelle
- Nach einem Austausch der Lampe (durchgeführt von einem Servicetechniker)
- Wenn die Wellenlängenvalidierung fehlschlägt

## Vorbereitungen

Sowohl für die Wellenlängenvalidierung als auch für die Raman-Wellenlängenkalibrierung *müssen* folgende Bedingungen erfüllt sein:

- Das System ist äquilibriert und die Umgebungsbedingungen sind stabil.
- Es befinden sich keine Luftblasen in der Messzelle. Es wird empfohlen, das Wasser zu entgasen. Stellen Sie eine Flussrate von 0,5 mL/min ein und warten Sie ca. 15 Minuten, bevor Sie beginnen, bis die Basislinie ausreichend stabil ist.
- Das Wasser weist keine Verunreinigungen auf. Verwenden Sie nur Wasser in Fluoreszenz-Qualität und verwenden Sie ggf. einen Filter oder eine geeignete Säule vor dem Detektor.

## Um eine Wellenlängenkalibrierung durchzuführen

Sie können die Wellenlängenkalibrierung über Chromeleon durchführen. Die Wellenlängenkalibrierung kann einige Minuten in Anspruch nehmen. In dieser Zeit kann keine Datenaufnahme erfolgen.

1. Führen Sie das Kommando **CalibrateRaman** aus, um die Kalibrierung zu starten.

Folgende Meldung erscheint in Chromeleon:

### Make sure that

- the system is equilibrated and conditions are stable
- fluorescence-grade water is flowing through the cell
- the water is degassed (no air bubbles) and free of particles
- Bestätigen Sie die Meldung mit OK, wenn Sie sicher sind, dass die oben genannten Bedingungen erfüllt sind.
   Die Kalibrierung kann bis zu fünf Minuten in Anspruch nehmen. Die neuen Kalibrierwerte werden im Detektor gespeichert.
- 3. Führen Sie eine Wellenlängenvalidierung durch.

## Sehen Sie dazu auch

Durchführung einer Wellenlängenvalidierung (> Seite 139)

## 7.6 Durchführung einer Wellenlängenvalidierung

Sie können die Wellenlängenkalibrierung mit Wasser (Raman-Messung) überprüfen. Schlägt die Validierung fehl, d.h. wurden deutliche Abweichungen in den Kalibrierwerten festgestellt, können Sie eine Wellenlängenkalibrierung mit Wasser (Raman-Messung) durchführen. Das Lichtspektrum der Xenon-Blitzlampe wird zur Kalibrierung der Anregungswellenlänge verwendet.

Wann

- Im Anschluss an die (Re)Installation des Detektors
- Im Anschluss an die Wellenlängenkalibrierung

#### Vorbereitungen

Sowohl für die Wellenlängenvalidierung als auch für die Raman-Wellenlängenkalibrierung *müssen* folgende Bedingungen erfüllt sein:

- Das System ist äquilibriert und die Umgebungsbedingungen sind stabil.
- Es befinden sich keine Luftblasen in der Messzelle. Es wird empfohlen, das Wasser zu entgasen. Stellen Sie eine Flussrate von 0,5 mL/min ein und warten Sie ca. 15 Minuten, bevor Sie beginnen, bis die Basislinie ausreichend stabil ist.
- Das Wasser weist keine Verunreinigungen auf. Verwenden Sie nur Wasser in Fluoreszenz-Qualität und verwenden Sie ggf. einen Filter oder eine geeignete Säule vor dem Detektor.

### So führen Sie eine Wellenlängenvalidierung durch

Sie können die Wellenlängenvalidierung über Chromeleon durchführen. Die Wellenlängenvalidierung kann einige Minuten in Anspruch nehmen. In dieser Zeit kann keine Datenaufnahme erfolgen.

1. Führen Sie den Befehl **ValidateRaman** aus, um die Validierung zu starten.

Folgende Meldung erscheint in Chromeleon:

Make sure that

- the system is equilibrated and conditions are stable
- fluorescence-grade water is flowing through the cell
- the water is degassed (no air bubbles) and free of particles

- Bestätigen Sie die Meldung mit OK, wenn Sie sicher sind, dass die oben genannten Bedingungen erfüllt sind.
   Die Wellenlängenvalidierung kann bis zu fünf Minuten in Anspruch nehmen. Im Chromeleon Audit Trail wird das Ergebnis ("passed"/"failed") angezeigt.
  - Passed: Die Wellenlängengenauigkeit ist innerhalb der Spezifikation. Die genaue gemessene Wellenlänge können Sie unter RamanValidationWL ablesen. Der Idealwert der Messung liegt bei 397 nm.
  - Failed: Die Kalibrierung des Detektors ist nicht ausreichend.
     Führen Sie eine Wellenlängenkalibrierung durch. Wenn die Wellenlängenvalidierung auch nach einer erneuten Kalibrierung fehlschlägt, prüfen Sie die unten aufgeführten Ursachen und Abhilfemaßnahmen.

Mögliche Ursache	Abhilfe
Messzelle falsch eingebaut	Prüfen Sie, ob die Messzelle korrekt sitzt und ob die Schrauben fingerfest angezogen sind.
Im System herrschen nicht die erforderlichen Bedingungen.	Beachten Sie die Bedingungen für die Wellenlängenvalidierung und -kalibrierung. Wird der Vorgang wiederholt abgebrochen, wenden Sie sich an den Kundendienst.
Messzelle kontaminiert	Reinigen Sie die Messzelle.

### Sehen Sie dazu auch

- Durchführung einer Wellenlängen-Kalibrierung () Seite 137)
- Reinigen der Messzelle (> Seite 144)

## 7.7 Messzelle

In diesem Abschnitt finden Sie Informationen zum Reinigen und Tausch von Messzellen. Zum Aus- und Einbau der Messzelle wird kein Werkzeug benötigt.

## 7.7.1 Hinweise zum Gebrauch von Messzellen

## ACHTUNG

Die Messzellen reagieren äußerst empfindlich auf Verschmutzung und Staub. Beachten Sie folgende Hinweise beim Gebrauch von Messzellen:

- Berühren Sie nie den optischen Block der Messzelle oder die empfindliche Elektronik auf der Rückseite.
- Die optischen Ports an der Messzelle sind empfindlich gegen Verunreinigungen und Kratzern. Berühren Sie niemals die optischen Ports der Messzelle und tauchen Sie diese niemals unter.
- Um Schäden an den optischen Ports der Messzelle zu vermeiden, gehen Sie vorsichtig vor, wenn Sie die Messzelle in den Messzellenschacht des Geräts einschieben.
- Auf der Rückseite der Messzelle befinden sich die empfindlichen Kontaktpunkte für den Identifikationschip. Berühren Sie niemals die Kontaktpunkte. Vermeiden Sie Schäden an der Elektronik des ID-Chips.
- Schmutzpartikel, Staub und Ablagerungen können zu einer Verunreinigung und Verstopfung der Messzelle führen. Wenn Kapillaren vom Einlass und Auslass der Messzelle entfernt werden, schließen Sie die Ports mit den Verschlussstopfen, um ein Verstopfen der Messzelle durch Schmutzpartikel bei der nächsten Anwendung zu vermeiden. Verwenden Sie die Stopfen, die bei Auslieferung an der Zelle installiert waren.
- Verwenden Sie die vorgesehene Verpackung f
  ür die Messzelle, wenn Sie sie aufbewahren oder transportieren.

## 7.7.2 Ausbauen der Messzelle

Erforderliche Teile

- Messzellen-Verpackung
- Abdeckung für den Messzellenschacht am Gerät

### Vorbereitungen

- 1. Spülen Sie die Messzelle beispielsweise mit Isopropanol, um sie von Lösungsmitteln frei zu spülen.
- 2. Schalten Sie den Pumpenfluss ab.

Gehen Sie wie folgt vor



## VORSICHT—Heiße Oberfläche

Die Messzelle kann sich erhitzen. Das Berühren einer heißen Messzelle kann zu Verbrennungen führen.

- Bevor Sie die Messzelle herausnehmen, berühren Sie die Messzelle kurz vorsichtig, um herauszufinden, ob sie heiß ist.
- Wenn die Messzelle heiß ist, lassen Sie die Messzelle abkühlen, bevor Sie die Messzelle herausnehmen. Stellen Sie sicher, dass die Temperatursteuerung für die Messzelle ausgeschaltet ist.
- 1. Entfernen Sie die Kapillaren vom Einlass und Auslass der Messzelle.
- 2. Verschließen Sie den Einlass und Auslass der Messzelle mit den Verschlussstopfen, die bei Auslieferung an der Messzelle installiert waren.
- 3. Lösen Sie die beiden Befestigungsschrauben der Messzelle. Die Schrauben sind unverlierbar mit der Messzelle verbunden und müssen nicht entfernt werden.
- 4. Entfernen Sie die Messzelle vom Messzellenschacht.

5. Montieren Sie die Abdeckung am Messzellenschacht des Detektors. Ziehen Sie die Schrauben fingerfest an (kein Werkzeug benötigt).



Abbildung 36: Einsetzen der Messzellenabdeckung

Nr.	Beschreibung
1	Befestigungsschrauben der Messzelle
2	Messzellen-Abdeckung

## ACHTUNG

Der Messzellenschacht am Gerät ist empfindlich gegen Staub und Schmutzpartikel.

Wenn keine Messzelle im Gerät eingebaut ist, verschließen Sie den Messzellenschacht mit der dafür vorgesehenen Abdeckung.

6. Lagern Sie die Messzelle zur Aufbewahrung und zum Versand in der Original-Verpackung.

## 7.7.3 Reinigen der Messzelle

Wann

Wenn Sie Ablagerungen von Eluenten- oder Probenbestandteilen an den Fenstern der Messzelle vermuten.

Erforderliche Teile

- Spül- und Injektionskit für Messzellen (optional)
- Wasser in HPLC-Qualität
- 0,1 molare Salpetersäure



## WARNUNG—Gesundheitsrisiken

Der Umgang mit Lösemitteln kann Gesundheits- und Sicherheitsrisiken darstellen.

Tragen Sie persönliche Schutzausrüstung je nach Gefahr und folgen Sie der Guten Laborpraxis. Informieren Sie sich im Sicherheitsdatenblatt des Herstellers.

## Gehen Sie wie folgt vor

Sie können die folgende Prozedur mit Hilfe des optional erhältlichen Spül- und Injektionskits durchführen.

- 1. Spülen Sie die Messzelle mit Wasser in HPLC-Qualität.
- 2. Füllen Sie die Messzelle mit Hilfe des optional erhältlichen Spül- und Injektionskits mit 0,1 molarer Salpetersäure.
- 3. Spülen Sie die Messzelle so lange mit Wasser in HPLC-Qualität, bis das Lösungsmittel aus der Messzelle wieder neutral (pH 7) ist.
- 4. Tauschen Sie die Messzelle aus, wenn sich das Problem nicht durch Reinigen beheben lässt.
### 7.7.4 Einbauen der Messzelle

Erforderliche Teile

Messzelle

Gehen Sie wie folgt vor

1. Setzen Sie die Messzelle gerade in den Messzellenschacht ein.



Abbildung 37: Einsetzen der Messzelle

- 2. Ziehen Sie die Befestigungsschrauben der Messzelle handfest an.
- Schließen Sie die Kapillaren an die Messzelle an. Folgen Sie den Anweisungen unter Anschließen der Flussverbindungen (▶ Seite 64). Stellen Sie den Fluss an und überprüfen Sie alle Verbindungen auf deren Dichtigkeit.
- 4. Schließen Sie die Gerätetüren, um die Erkennung der Messzelle zu ermöglichen. Alternativ können Sie ebenso die **INIT**-Taste auf der Tastatur drücken.
- 5. Führen Sie eine Wellenlängenkalibrierung durch.

#### Sehen Sie dazu auch

Durchführung einer Wellenlängen-Kalibrierung () Seite 137)

# 7.8 Tauschen der Sicherungen

#### Wann

Durchgebrannte Sicherungen

#### Erforderliche Teile

Sicherungen (2 Sicherungen, 3.15 AT, 250 V AC, träge, 5 x 20 mm) aus dem Sicherungskit

#### Erforderliche Werkzeuge

Schlitzschraubendreher, jede Größe zwischen 3,3 mm und 5,5 mm ist geeignet

#### Vorbereitungen



#### WARNUNG—Stromschlag

Im Gerät treten hohe Spannungen auf, die zu einem Stromschlag führen oder Schäden am Gerät verursachen können.

- Schalten Sie das Gerät über den Hauptnetzschalter aus. Trennen Sie das Netzkabel von der Stromversorgung und vom Gerät.
- Verwenden Sie nur Sicherungen des Typs und der Stromstärke, die von Thermo Fisher Scientific für das Gerät spezifiziert sind.
- Verwenden Sie keine reparierten Sicherungen und schließen Sie den Sicherungshalter nicht kurz.

#### Gehen Sie wie folgt vor

Der Sicherungshalter befindet sich neben dem Hauptnetzschalter.



Abbildung 38: Sicherungshalter

Nr.	Beschreibung
1	Hauptnetzschalter (Ein/Aus)
2	Sicherungshalter
3	Netzbuchse

- 1. Entfernen Sie den Sicherungshalter mit dem Schraubendreher.
- 2. Tauschen Sie die beiden Sicherungen gegen Sicherungen des spezifizierten Typs und der spezifizierten Stromstärke aus. Tauschen Sie immer *beide* Sicherungen aus.
- 3. Setzen Sie den Sicherungshalter wieder ein.
- 4. Schließen Sie das Netzkabel wieder an die Stromversorgung und an das Gerät an.
- 5. Schalten Sie das Gerät über den Hauptnetzschalter ein.

# 7.9 Aktualisieren der Gerätefirmware

#### Wann

Eine Aktualisierung der Gerätefirmware kann erforderlich sein, zum Beispiel, wenn es eine neue Firmware-Version gibt, die neue Funktionen enthält oder Probleme einer früheren Version behebt.

#### Erforderliche Utensilien

Firmware-Version/Chromeleon-Version, wie erforderlich

**TIPP** Wenn eine neue Firmware-Version verfügbar ist, wird diese zusammen mit der nächsten Chromeleon-Version ausgeliefert. Die neue Firmware wird *nicht* automatisch auf das Gerät übertragen, wenn Sie die Chromeleon-Version installieren.

#### Vorbereitungen

- 1. Lesen Sie die Release-Notes, die mit der Firmware und/oder Chromeleon-Version mitgeliefert werden.
- 2. Verbinden Sie das Gerät in der Software Chromeleon.
- 3. Halten Sie alle Operationen auf der Anlage (englisch: Instrument) an, die das Gerät enthält.
- Warten Sie, bis die Anlage nicht mehr in Verwendung ist (Modus: 'idle').

#### Gehen Sie wie folgt vor

- 1. Starten Sie das Programm Instrument Configuration Manager.
- Aktualisieren Sie die Firmware über die Registerkarte General im Konfigurationsdialog für das Gerät. Einzelheiten hierzu finden Sie in der Chromeleon-Hilfe.
   Die Aktualisierung der Firmware kann einige Minuten dauern.

#### ACHTUNG

Das Aufspielen einer früheren Firmware-Version oder eine nicht vollständig ausgeführte Aktualisierung der Firmware kann zu Funktionsverlusten oder Fehlfunktionen des Geräts führen.

- Unterbrechen Sie während des Downloads keinesfalls die Kommunikation zwischen der Software Chromeleon und dem Gerät.
- Zu Beginn der Aktualisierung informiert Sie eine Meldung über die aktuell im Gerät installierte Firmware-Version und die Version, die von der Software Chromeleon auf das Gerät übertragen wird. Brechen Sie die Installation ab, wenn die im Gerät installierte Firmware neuer ist als die Version in der Software Chromeleon.
- 3. Überprüfen Sie im Audit Trail des Programms Instrument Configuration Manager, ob die Firmware-Aktualisierung erfolgreich war oder fehlgeschlagen ist.
  - Wenn die Aktualisierung der Firmware fehlgeschlagen ist, schalten Sie das Gerät aus und wieder ein und wiederholen Sie die Aktualisierung.
  - Wenn die Aktualisierung der Firmware wiederholt fehlschlägt, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst.
- Nach erfolgreicher Aktualisierung der Firmware, kann eine Requalifizierung des Geräts erforderlich sein. Empfehlungen finden Sie in den Release Notes.

# 7.10 Tauschen der Türen

#### Wann

Beschädigung der Tür

**TIPP** Es ist nicht erforderlich, die Türen für die Wartungsarbeiten zu entfernen. Sollte dies aus einem bestimmten Grund oder für eine bestimmte Arbeit erforderlich sein, folgen Sie den entsprechenden Schritten in diesem Abschnitt.

#### Erforderliche Teile

Ersatztür

Gehen Sie wie folgt vor

#### ACHTUNG

Um die Türangeln nicht zu beschädigen, gehen Sie bei den folgenden Schritten vorsichtig vor und wenden Sie keine Gewalt an.

1. Zum Entfernen einer Tür drücken Sie die Tür beim Öffnen nach oben.

Öffnen Sie die Tür bis zu einer Position, in der die beiden Türangeln am Gehäuse mit den Nuten in der Tür ausgerichtet sind. Sie können die Tür nur entfernen, wenn sich die Angeln in den Nuten befinden.



Abbildung 39: Aushängen einer Tür

Nr.	Beschreibung
1	Türangel am Gehäuse
2	Aufnahmenut der Tür

2. Kippen Sie die Tür leicht nach außen, weg vom Gehäuse, und entfernen Sie die Tür.

- 3. Zum Einbau der Tür richten Sie die Tür mit den Türangeln am Gehäuse aus. Achten Sie darauf, keine Schläuche oder Kapillaren zwischen Tür und Gehäuse einzuklemmen.
- 4. Setzen Sie die Türangeln in die Nut ein, indem Sie die Tür nach oben drücken und leicht drehen.
- Drücken Sie die Tür nach unten, damit sie einrastet. Sie können die Tür nur schließen, wenn sie korrekt eingebaut ist.

# 7.11 Transportieren und Versenden des Gerätes

Wenn Sie das Gerät an einen anderen Ort transportieren möchten oder versenden müssen, bereiten Sie das Gerät zuerst für den Transport vor und transportieren oder versenden Sie das Gerät dann wie erforderlich. Folgen Sie den Anweisungen in diesem Abschnitt.

Beachten Sie folgende Sicherheitshinweise:



#### VORSICHT—Schweres und unhandliches Gerät

Das Gerät ist zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden am Gerät zu vermeiden, beachten Sie folgende Hinweise:

- Um das Gerät zu greifen und zu transportieren, sind mindestens zwei Personen erforderlich, zum Beispiel, um das Gerät anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn das Gerät in den Systemturm gehoben oder aus dem Systemturm entfernt wird.
- Verwenden Sie f
  ür den Transport des Ger
  äts die Tragegriffe, die mitgeliefert wurden. Transportieren Sie oder heben Sie das Ger
  ät niemals an den Ger
  ätet
  üren. Dies f
  ührt zu Besch
  ädigungen an den T
  üren oder am Ger
  ät.

#### Gehen Sie wie folgt vor

- 1. Bereiten Sie das Gerät für den Transport vor. Siehe Vorbereiten des Gerätes für den Transport (▶ Seite 152).
- 2. Dieser Schritt ist wie folgt abhängig:
  - Um das Gerät an einen anderen Standort zu transportieren, folgen Sie den Anweisungen in Transportieren des Geräts an einen anderen Standort (> Seite 153).
  - ◆ Um das Gerät zu versenden, folgen Sie den Anweisungen in Versenden des Gerätes (▶ Seite 154).

#### 7.11.1 Vorbereiten des Gerätes für den Transport

Bereiten Sie das Gerät wie folgt für den Transport vor:

 Führen Sie die Maßnahmen zur langfristigen Außerbetriebnahme des Geräts durch (siehe Langfristige Außerbetriebnahme (▶ Seite 122)).

- 2. Schalten Sie das Gerät über den Hauptnetzschalter aus und ziehen Sie den Netzstecker.
- 3. Entfernen Sie alle Kabel und Flussverbindungen zu anderen Geräten.
- Bauen Sie die Messzelle aus Detektor aus und bewahren Sie diese in ihrer Verpackung auf. Achten Sie darauf, dass Sie die Abdeckung am Messzellenschacht montieren. Messzellen müssen in ihrer original Messzellen-Verpackung versendet werden.

#### ACHTUNG

Der Messzellenschacht am Gerät ist empfindlich gegen Staub und Schmutzpartikel.

Wenn keine Messzelle im Gerät eingebaut ist, verschließen Sie den Messzellenschacht mit der dafür vorgesehenen Abdeckung.

 Entfernen Sie, je nach Erfordernis, das Gerät oder den Moduleinschub aus dem Systemturm: Montieren Sie die Tragegriffe und entfernen Sie das Gerät aus dem Vanquish-System. Informationen zum Abbau des Systemturms finden Sie im Abschnitt Transportieren oder Versenden des Systems in der Vanquish-System-Betriebsanleitung.

-oder-

Entfernen Sie den Moduleinschub aus dem Gerätegehäuse im Systemturm (siehe Entfernen des Moduleinschubs (▶ Seite 156)).

#### Sehen Sie dazu auch

Ausbauen der Messzelle (▶ Seite 142)

#### 7.11.2 Transportieren des Geräts an einen anderen Standort

#### Vorbereitungen

Bereiten Sie das Gerät für den Transport vor. Siehe Vorbereiten des Gerätes für den Transport (▶ Seite 152).

#### Gehen Sie wie folgt vor

- 1. Beachten Sie die Hinweise zum sicheren Heben und Bewegen des Gerätes.
- 2. Transportieren Sie das Gerät an den neuen Standort.
- 3. Installieren und schließen Sie das Gerät im Systemturm an. Folgen Sie der Anweisung zum Aufbau des Systemturms in der *Vanquish-System-Betriebsanleitung*.

- 4. Schließen Sie das Gerät an:
  - a) Verbinden Sie das Gerät und stellen Sie die fluidischen Verbindungen her (siehe Installation (▶ Seite 47)).
  - b) Bereiten Sie das Gerät für den Betrieb vor (siehe Vorbereiten des Gerätes für den Betrieb (▶ Seite 98)).
- 5. Bevor Sie mit einer Analyse beginnen, warten Sie, bis sich das Gerät äquilibriert hat, und vergewissern Sie sich, dass es für den Betrieb bereit ist.

### 7.11.3 Versenden des Gerätes

#### Vorbereitungen

Bereiten Sie das Gerät für den Transport vor. Siehe Vorbereiten des Gerätes für den Transport (▶ Seite 152).

#### Gehen Sie wie folgt vor

- Folgen Sie den Anweisungen zum Auspacken in dieser Anleitung in umgekehrter Reihenfolge.
   Verwenden Sie ausschließlich die Original-Verpackungsmaterialien und Originalverpackung. Ist die Originalverpackung nicht mehr verfügbar, können Sie geeignete Geräteverpackungen über die Thermo Fisher Scientific-Vertriebsorganisation bestellen.
- Wenden Sie sich an den Thermo Fisher Scientific-Kundendienst vor Ort, wenn Sie das Gerät für eine Reparatur an Thermo Fisher Scientific zurückschicken müssen.



#### VORSICHT—Gefahr durch Kontamination

Das Gerät wurde möglichweise während des Betriebs mit gefährlichen Substanzen kontaminiert, welche die Gesundheit des Servicepersonals gefährden können.

- Dekontaminieren Sie alle Teile des Geräts, die Sie zur Reparatur zurückschicken möchten.
- Füllen Sie das Gesundheits- und Sicherheitsformular ("Health and Safety Form") aus. Thermo Fisher Scientific nimmt keine Geräte zur Reparatur an, wenn das Gesundheits- und Sicherheitsformular fehlt, unvollständig ausgefüllt ist, oder nicht unterschrieben ist.

#### Wiederaufnehmen des Gerätebetriebs nach einem Versand

Um das Gerät nach einem Versand wieder in Betrieb zu nehmen, gehen Sie wie folgt vor:

- 1. Folgen Sie den Anweisungen zum Auspacken in dieser Betriebsanleitung.
- 2. Installieren und schließen Sie das Gerät im Systemturm an. Folgen Sie der Anweisung zum Aufbau des Systemturms in der *Vanquish-System-Betriebsanleitung*.
- 3. Schließen Sie das Gerät an:
  - a) Verbinden Sie das Gerät und stellen Sie die fluidischen Verbindungen her (siehe Installation (▶ Seite 47)).
  - b) Bereiten Sie das Gerät für die Erstinbetriebnahme vor (siehe Vorbereiten des Gerätes für den Betrieb (▶ Seite 98)).
- 4. Bevor Sie mit einer Analyse beginnen, warten Sie, bis sich das Gerät äquilibriert hat, und vergewissern Sie sich, dass es für den Betrieb bereit ist.

# 7.12 Tauschen des Moduleinschubs

### 7.12.1 Entfernen des Moduleinschubs



#### VORSICHT—Schweres und unhandliches Gerät

Das Gerät ist zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden am Gerät zu vermeiden, beachten Sie folgende Hinweise:

- Um das Gerät zu greifen und zu transportieren, sind mindestens zwei Personen erforderlich, zum Beispiel, um das Gerät anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn das Gerät in den Systemturm gehoben oder aus dem Systemturm entfernt wird.

Erforderliche Werkzeuge

Schraubendreher, Torx T20

Vorbereitungen

1. Bereiten Sie das Gerät für den Transport vor. Siehe Transportieren und Versenden des Gerätes (▶ Seite 152).

#### Gehen Sie wie folgt vor

1. Lösen Sie die vier Unverlierbarkeitsschrauben vorne links und vorne rechts am Gerät.



Abbildung 40: Unverlierbarkeitsschrauben am Moduleinschub

 Drücken Sie alle Schläuche und Kapillaren, die sich im Schlauchkanal der Vanquish-Systemmodule befinden, in den Schlauchkanal hinein. Andernfalls können Sie den Moduleinschub im nächsten Schritt nicht richtig aus dem Gehäuse entfernen.  Greifen Sie den Moduleinschub am Panel unterhalb des Messzellenschachts oder an der Leckage-Wanne und ziehen Sie das Modul circa 10 cm aus dem Gehäuse heraus.



Abbildung 41: Herausziehen des Moduleinschubs

#### ACHTUNG

Der Moduleinschub kann herunterfallen, wenn Sie ihn zu weit aus dem Gehäuse herausziehen.

Ziehen Sie den Moduleinschub nur so weit hinaus, dass Sie ihn an beiden Seiten von unten greifen können.

- 4. Entfernen Sie den Moduleinschub aus dem Gehäuse. Für die folgenden Schritte sind mindestens zwei Personen erforderlich:
  - a) Greifen Sie den Moduleinschub auf beiden Seiten von unten.
  - b) Ziehen Sie den Moduleinschub von den Schienen nach vorn heraus.
  - c) Stellen Sie den Moduleinschub auf eine saubere und stabile Oberfläche.

### 7.12.2 Zurückschicken des Moduleinschubs

#### Vorbereitung

*Falls noch nicht erfolgt:* Entfernen Sie den Moduleinschub aus dem Gehäuse. Siehe Entfernen des Moduleinschubs (▶ Seite 156).

#### Gehen Sie wie folgt vor

- Bauen Sie den Moduleinschub in das Gehäuse der Ersatzteilgeräts ein. Folgen Sie den Schritten für den Einbau des Moduleinschubs in das Gehäuse und zum Anziehen der Unverlierbarkeitsschrauben im Installieren des Moduleinschubs (> Seite 159).
- 2. Folgen Sie den Anweisungen in Versenden des Gerätes (> Seite 154).

#### ACHTUNG

Wird der Moduleinschub nicht ordnungsgemäß verschickt, führt dies zu Schäden am Gerät.

Versenden Sie den Moduleinschub immer entsprechend der Beschreibung in dieser Betriebsanleitung.



#### VORSICHT—Gefahr durch Kontamination

Das Gerät wurde möglichweise während des Betriebs mit gefährlichen Substanzen kontaminiert, welche die Gesundheit des Servicepersonals gefährden können.

- Dekontaminieren Sie alle Teile des Geräts, die Sie zur Reparatur zurückschicken möchten.
- Füllen Sie das Gesundheits- und Sicherheitsformular ("Health and Safety Form") aus. Thermo Fisher Scientific nimmt keine Geräte zur Reparatur an, wenn das Gesundheits- und Sicherheitsformular fehlt, unvollständig ausgefüllt ist, oder nicht unterschrieben ist.

### 7.12.3 Installieren des Moduleinschubs



#### VORSICHT—Schweres und unhandliches Gerät

Das Gerät ist zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden am Gerät zu vermeiden, beachten Sie folgende Hinweise:

- Um das Gerät zu greifen und zu transportieren, sind mindestens zwei Personen erforderlich, zum Beispiel, um das Gerät anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn das Gerät in den Systemturm gehoben oder aus dem Systemturm entfernt wird.

#### Erforderliche Werkzeuge

Schraubendreher, Torx T20

#### Vorbereitungen

- Packen Sie das Ersatzgerät aus. Entfernen Sie den Moduleinschub aus dem Gehäuse wie in Entfernen des Moduleinschubs (>> Seite 156) beschrieben.
- Vergewissern Sie sich, dass das Gerätegehäuse, in das der Moduleinschub installiert werden soll, sauber ist. Reinigen Sie gegebenenfalls die Gehäuseoberflächen von innen und von außen. Siehe Reinigen oder Dekontaminieren des Gerätes (> Seite 133).
- Wenn Sie den Moduleinschub in ein Gehäuse im Systemturm installieren, kontrollieren Sie, dass das Gehäuse ordnungsgemäß im Turm platziert ist.

#### Gehen Sie wie folgt vor

- Drücken Sie alle Schläuche und Kapillaren, die sich im Schlauchkanal der Vanquish-Systemmodule befinden, in den Schlauchkanal hinein. Andernfalls können Sie den Moduleinschub im nächsten Schritt nicht richtig in das Gehäuse einsetzen.
- Prüfen Sie am Moduleinschub, ob Sie die Unverlierbarkeitsschrauben mit den Fingern hin und her bewegen können. Wenn Sie dies nicht können, schrauben Sie sie soweit ein, bis Sie sie es können.

- 3. Setzen Sie den Moduleinschub in das Gehäuse ein. Für die folgenden Schritte sind mindestens zwei Personen erforderlich:
  - a) Greifen Sie den Moduleinschub auf beiden Seiten von unten.
  - b) Heben Sie den Moduleinschub auf Gehäusehöhe an.
  - c) Platzieren Sie den Moduleinschub so im Gehäuse, dass sich das Modul circa 25 cm weit im Gehäuse befindet.
  - d) Schieben Sie den Moduleinschub auf die Schienen und in das Gehäuse, bis sich das Modul vollständig im Gehäuse befindet.



Abbildung 42: Einsetzen des Moduleinschubs

- 4. Damit Sie den Moduleinschubs leichter im Gehäuse befestigen können, ziehen Sie den Moduleinschub einige Millimeter heraus (circa 2 bis 5 mm).
- 5. Bewegen Sie jede Unverlierbarkeitsschraube mit dem Schraubendreher hin und her und drücken Sie sie dabei nach innen, bis die Schraube in das Gewinde gleitet.
- 6. Schieben Sie den Moduleinschub bis zum Anschlag in das Gehäuse zurück.
- 7. Ziehen Sie die vier Unverlierbarkeitsschrauben des Moduleinschubs gleichmäßig und schrittweise handfest an.

#### ACHTUNG

- Vergewissern Sie sich, dass die Schrauben fest angezogen sind.
   Ziehen Sie den Moduleinschub an der Leckage-Wanne nach vorn und kontrollieren Sie, ob sich die Schrauben bewegen. Bewegen Sie sich nicht, ist der Moduleinschub ordnungsgemäß installiert.
- Wenn sich die Schrauben bewegen, ziehen Sie diese nach. Wird ein Drehmomentschraubendreher verwendet, ist das empfohlene Drehmoment 1,2 Nm.
- 8. Schließen Sie das Gerät an und nehmen Sie es wieder in Betrieb.

### 7.12.4 Anschließen des Moduleinschubs

Wenn Sie den Moduleinschub wieder in das Gehäuse eingesetzt haben, schließen Sie das Gerät an und nehmen Sie es wieder in Betrieb.

#### Gehen Sie wie folgt vor

- 1. Schließen Sie den Moduleinschub an:
  - a) Verbinden Sie den Moduleinschub und stellen Sie die fluidischen Verbindungen her (siehe Installation (▶ Seite 47)).
- 2. Bereiten Sie *alle anderen* Module des Vanquish-Systems für den Betrieb vor und nehmen Sie sie in Betrieb. Informieren Sie sich in den *Betriebsanleitungen* der anderen Module.
- 3. Bevor Sie mit einer Analyse beginnen, warten Sie, bis sich das Chromatographie-System äquilibriert hat, und vergewissern Sie sich, dass es für den Betrieb bereit ist.
- 4. Führen Sie in der Software Chromeleon die gerätespezifischen **Performance Qualification** (PQ) Tests durch.
- 5. Wenn der Test abgeschlossen ist, aktualisieren Sie den Parameter **QualificationDone**.

# 8 Fehlersuche

In diesem Kapitel finden Sie einen Leitfaden zur Fehlersuche für den Betrieb des Gerätes.

# 8.1 Allgemeine Informationen zur Fehlersuche

Bei der Erkennung und Behebung von Störungen, die beim Betrieb des Gerätes auftreten können, helfen Ihnen nachfolgende Funktionen.

**TIPP** Informationen zu Störungen, die beim Betrieb eines Vanquish-Systems auftreten können, finden Sie in der *Vanquish-System-Betriebsanleitung*.

Wenn Sie eine Störung mithilfe der Beschreibungen in diesem Kapitel nicht selbst beheben können, oder Sie auf Probleme stoßen, die hier nicht beschrieben sind, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst, um Unterstützung zu erhalten. Kontaktinformationen finden Sie am Anfang dieser Betriebsanleitung.

Für die Kommunikation mit Thermo Fisher Scientific benötigen Sie zur leichteren Identifizierung Ihres Gerätes die Seriennummer und den technischen Namen.

#### Statusanzeigen

Die LED-Statusleiste an der Vorderseite des Gerätes sowie die **STATUS**-LED auf der Tastatur im Gerät ermöglichen einen schnellen, visuellen Überblick über den Betriebszustand des Geräts. Wenn die Geräte-Firmware ein Problem erkennt, leuchten die Statusanzeigen rot. Das Problem wird an die Software Chromeleon weitergereicht und es erscheint eine Meldung im Audit Trail.

#### Alarme

Undichtigkeiten sind ein potentielles Sicherheitsrisiko. Wenn der Leaksensor auf eine Undichtigkeit anspricht, ertönt daher zusätzlich zu der Meldung im Chromeleon Audit Trail ein akustisches Signal und die Statusanzeige leuchtet rot. Finden und beseitigen Sie die Ursache wie in diesem Handbuch beschrieben.

Meldungen im Chromeleon Audit Trail

Wenn die Geräte-Firmware ein Problem erkennt, wird dieses an die Software Chromeleon weitergereicht.

Die Software Chromeleon protokolliert Informationen über alle Ereignisse, welche die Gerätesteuerung betreffen, tagesweise in einem Audit Trail (Protokoll). Der Audit Trail wird nach dem Datum benannt, im Datumsformat yyyymmdd (jjjjmmtt). Der Audit Trail für den 15. Mai 2019 ist also 20190515 benannt. Die Anlagenprotokolle (Instrument Audit Trails) befinden sich auf dem ePanel Set (ePanel "Audit"). Zusätzlich stehen Audit Trails für jede Anlage in der Datenansicht der Console im Ordner der jeweiligen Anlage zur Verfügung.

Den Meldungen im Chromeleon Audit Trail ist ein Symbol vorangestellt. Das Symbol zeigt den Schweregrad des Problems an (siehe *Chromeleon-Hilfe*). Mögliche Fehlerursachen und empfohlene Abhilfemaßnahmen finden Sie in Meldungen (▶ Seite 166).

# 8.2 Meldungen

Die Tabelle führt die am häufigsten beobachteten Meldungen für das Gerät auf und leistet Hilfe bei der Fehlersuche und Fehlerbehebung.

Jede Meldung besteht aus einer Code-Nummer und einem Text. Die Code-Nummer ist das eindeutige Identifizierungszeichen, während sich die Formulierung des Textes verändern kann. Beachten Sie Folgendes:

- Zum leichteren Auffinden einer Meldung, erscheinen die Meldungen in der Tabelle nach Code sortiert.
- Wenn Sie einen Code nicht finden können, pr
  üfen Sie den Text der Meldung. Die beiden Meldungen "Unexpected module behavior" und "Module malfunction detected" können verschiedenen Codes zugewiesen sein. Sie finden eine Beschreibung am Anfang der Tabelle.

**TIPP** Wenn Sie die Störung nicht entsprechend der Anweisungen in dieser Anleitung beheben können oder wenn ein Code oder eine Meldung erscheint, die nicht in der Tabelle aufgeführt ist, notieren Sie den Code und den Wortlaut der Meldung und wenden Sie sich an uns. Siehe dazu *Kontaktinformationen* am Anfang dieser Betriebsanleitung.

Meldung und Code	Beschreibung und Abhilfe
Unexpected module behavior. Code xx	Unerwartetes Modulverhalten. xx = Zwei- bis vierstellige Code-Nummer. Wenn die Meldung erscheint, notieren Sie den Code der Meldung und schalten Sie das Gerät aus. Warten Sie 5 Sekunden und schalten Sie das Gerät wieder ein. <b>TIPP</b> Wenn die Meldung mit <i>Code 103</i> erscheint, sind weitere Abhilfemaßnahmen möglich. Siehe weiter unten in dieser Tabelle. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Module malfunction detected. Code xx	Fehlfunktion des Moduls erkannt. xx = Zwei- bis vierstellige Code-Nummer. Wenn die Meldung erscheint, notieren Sie den Code der Meldung. Schalten Sie das Gerät aus und wenden Sie sich an den Kundendienst.
Code 33 Leak detected – eliminate within approx. xx seconds.	Undichtigkeit erkannt - beheben Sie diese innerhalb von xx Sekunden. xx = Zeit in Sekunden, um die Undichtigkeit zu beheben Finden und beheben Sie die Ursache für die Undichtigkeit (siehe Beheben von Undichtigkeiten (▶ Seite 171)).
Code 34 Leak detected.	Undichtigkeit erkannt. Finden und beheben Sie die Ursache für die Undichtigkeit (siehe Beheben von Undichtigkeiten (▶ Seite 171)).

Meldung und Code	Beschreibung und Abhilfe
Code 36 Download failed.	Download fehlgeschlagen. Der Firmware-Download wurde nicht erfolgreich abgeschlossen. Prüfen Sie, ob Sie die richtige Firmware-Datei ausgewählt haben. Wiederholen Sie den Download.
Code 37 Download firmware mismatch.	Fehler bei Firmware-Download. Der Firmware-Download wurde nicht erfolgreich abgeschlossen. Prüfen Sie, ob Sie die richtige Firmware-Datei ausgewählt haben. Wiederholen Sie den Download.
Code 89 Liquid leak sensor missing or defective.	Leaksensor für Flüssigkeit fehlt oder ist defekt. Vergewissern Sie sich, dass der Leaksensor richtig installiert ist und die Kabelverbindung richtig angesteckt ist. Wenden Sie sich für Unterstützung an den Kundendienst von Thermo Fisher Scientific, wenn der Leaksensor defekt ist. Um das Gerät trotzdem zu betreiben, können Sie die Leaksensor- Funktionalität in der Software Chromeleon deaktivieren, indem Sie den Parameter <b>Leak Sensor Mode</b> auf <b>Disabled</b> setzen.
Code 90 Download firmware mismatch – invalid version.	Fehler bei Firmware-Download – ungültige Version. Sie haben versucht, eine frühere Firmware-Version aufzuspielen als aktuell im Gerät installiert ist. Das Aufspielen einer früheren Firmware-Version kann zu Funktionsverlusten oder Fehlfunktionen des Gerätes führen. Falls erforderlich, wiederholen Sie die Installation mit einer Firmware-Version, die neuer ist als die aktuell im Gerät installierte Version.
Code 103 Unexpected module behavior – limited features available.	Unerwartetes Modulverhalten - eingeschränkter Funktionsumfang. Die Firmware ist möglichweise defekt oder es wurde eine niedrigere Firmware-Version aufgespielt. Aktualisieren Sie die Firmware auf die neueste Version. Siehe Aktualisieren der Gerätefirmware (▶ Seite 148).
Code 118 USB Buffer Overflow.	<ul> <li>Überlauf des USB-Puffers.</li> <li>Es handelt sich um ein Software-Problem. Das Modul produziert Daten schneller als sie der Rechner, auf dem die Software Chromeleon läuft, lesen kann.</li> <li>1. Trennen Sie das Modul in der Software Chromeleon und verbinden Sie es erneut.</li> <li>2. Wird das Problem damit nicht gelöst, aktualisieren Sie die Firmware oder die Version der Software Chromeleon.</li> <li>3. Besteht das Problem weiterhin: Auch Software von Drittanbietern auf dem Rechner, zum Beispiel Virenscanner, können das Problem verursachen. Wenden Sie sich an die IT-Abteilung vor Ort.</li> </ul>
Code 120 The request timed out.	<ul> <li>Die Zeit für die Anforderung ist abgelaufen.</li> <li>Mit dem Modul kann keine Kommunikation aufgebaut werden. Das Modul hat nicht innerhalb der Zeit geantwortet.</li> <li>Für das Modul, für das die Meldung erscheint:</li> <li>1. Schalten Sie das Modul ein, wenn es noch nicht eingeschaltet ist.</li> <li>2. Überprüfen Sie, dass die System-Interlink-Kabel korrekt angeschlossen sind.</li> </ul>

Meldung und Code	Beschreibung und Abhilfe
Code 8013 No flow cell detected	Keine Messzelle erkannt Überprüfen Sie, ob eine Messzelle ordnungsgemäß installiert ist. Die Lampe wird automatisch ausgeschaltet, wenn keine Messzelle erkannt wurde.
Code 8032 Unexpected module behavior. Flow cell Code 8033 Unexpected module behavior. Flow cell	Unerwartetes Modulverhalten. Messzelle Wenn die Meldung "Flow cell" enthält: Überprüfen Sie, ob die Messzelle ordnungsgemäß installiert ist. Öffnen und schließen Sie die Türen. Schalten Sie den Detektor aus. Warten Sie 5 Sekunden und schalten Sie den Detektor wieder ein. Ersetzen Sie die Messzelle durch eine andere und versuchen Sie es erneut. Wenden Sie sich an den Kundendienst wenn die Meldung erneut erscheint
Code 8035 Signal overflow in <channel name=""></channel>	Signalüberlauf in <name des="" kanals=""> Das Referenzsignal ist zu gering oder das PMT-Signal ist zu groß. Prüfen Sie, ob die Messzelle kontaminiert ist. Reinigen Sie gegebenenfalls die Messzelle oder tauschen Sie diese aus. Überprüfen Sie, ob die Probenkonzentration zu hoch ist und stellen Sie sicher, dass die Anregungswellenlänge oberhalb von der UV- Cutoff-Wellenlänge des Eluenten eingestellt ist. Eventuell ist die Lampe zu alt und muss ausgetauscht werden.</name>
Code 8036 Reference signal too low in <channel name=""></channel>	Referenzsignal zu gering in <name des="" kanals=""> Das Referenzsignal ist zu gering (hohe Absorption). Prüfen Sie, ob die Messzelle kontaminiert ist. Reinigen Sie gegebenenfalls die Messzelle oder tauschen Sie diese aus. Überprüfen Sie, ob die Probenkonzentration zu hoch ist und stellen Sie sicher, dass die Anregungswellenlänge oberhalb von der UV-Cutoff-Wellenlänge des Eluenten eingestellt ist. Eventuell ist die Lampe zu alt und muss ausgetauscht werden.</name>
Code 8049 Command rejected - close front door first	Befehl abgelehnt - Schließen Sie zuerst die Gerätetür Es ist nicht möglich, die Datenaufnahme zu starten, wenn die Gerätetüren offen sind. Schließen Sie die Türen und versuchen Sie es erneut.
Code 8051 Flow cell detected. It is recommended executing the command CalibrateRaman	Messzelle erkannt. Es wird empfohlen, den Befehl CalibrateRaman auszuführen Es wurde eine neue Messzelle erkannt. Es wird empfohlen, eine Raman-Wellenlängenkalibrierung durchzuführen (siehe Durchführung einer Wellenlängen-Kalibrierung () Seite 137)), da ohne Kalibrierung die Wellenlängen bis zu einigen Nanometern von der tatsächlichen Wellenlänge abweichen können.
Code 8064 PMT x signal overflow in <channel name=""> Code 8065 PMT x signal overflow during scan</channel>	PMT x Signalüberlauf in <name des="" kanals=""> oder PMT x Signalüberlauf während des Scans Wobei x = PMT 1 oder 2 Signalüberlauf in PMT 1 oder 2 aufgetreten. Wiederholen Sie den Lauf oder Scan mit einer geringeren Sensitivity oder einer geringeren Probenkonzentration.</name>

Meldung und Code	Beschreibung und Abhilfe
Code 8071 PMT Sensitivity autorange for <channel name=""> -new Sensitivity is x</channel>	<ul> <li>PMT Empfindlichkeit Autorange für <name des="" kanals=""> - neue Empfindlichkeit ist x</name></li> <li>Wobei x = Die neue Empfindlichkeit nach der Anpassung.</li> <li>Die Empfindlichkeit musste automatisch reduziert werden. Wählen Sie eine geeignete Empfindlichkeit (siehe Sensitivity (Detektorempfindlichkeit) (▶ Seite 111)) und wiederholen Sie den Vorgang. x gibt die neue Empfindlichkeit nach der Anpassung an.</li> <li>Verwenden Sie diesen oder einen kleineren Wert in Ihrer Methode, wenn Sie ähnliche Konzentrationen erwarten.</li> </ul>
Code 8078 PMT Autorange for <channel name=""> not possible - already minimal Sensitivity</channel>	PMT Autorange für <name des="" kanals=""> nicht möglich, da bereits auf minimaler Empfindlichkeit Der Detektor konnte nicht automatisch die Empfindlichkeit reduzieren, da die Empfindlichkeit bereits auf ein Minimum eingestellt worden ist. Wiederholen Sie den Lauf oder Scan mit einer geringeren Probenkonzentration oder einem kleineren Injektionsvolumen.</name>
Code 8087 System Interlink error	System-Interlink Fehler Die System Interlink-Verbindung wurde unterbrochen. Prüfen Sie, dass alle System Interlink-Kabel am Vanquish-System korrekt angeschlossen sind.
Code 8093 Not initialized	Nicht initialisiert Der Detektor ist nicht initialisiert. Prüfen Sie, dass eine Messzelle korrekt installiert ist und dass sich die Eluentenzusammensetzung in der Messzelle nicht ändert und keine Luftblasen vorhanden sind. Öffnen und schließen Sie dann die Türen oder drücken Sie die <b>INIT</b> auf der Tastatur.
Code 8100 - Code 8109 Calibration of xx failed.	<ul> <li>Die Kalibrierung von xx ist fehlgeschlagen.</li> <li>Wobei x = Name der Komponente, für welche die Kalibrierung fehlgeschlagen ist.</li> <li><i>Mögliche Ursachen</i></li> <li>Die Raman-Wellenlängenkalibrierung ist fehlgeschlagen.</li> <li>Eventuell ist die Messzelle nicht korrekt eingebaut.</li> <li>Eventuell ist die Messzelle kontaminiert.</li> <li><i>Abhilfen</i></li> <li>Beachten Sie die Bedingungen für die Wellenlängenkalibrierung (siehe Durchführung einer Wellenlängen-Kalibrierung (▶ Seite 137)) und Wellenlängenvalidierung (siehe Durchführung einer Wellenlängenvalidierung (siehe 139)).</li> <li>Prüfen Sie, dass die Messzelle korrekt eingebaut ist und dass die Schrauben fingerfest angezogen sind.</li> <li>Reinigen Sie die Messzelle (siehe Reinigen der Messzelle (▶ Seite 144)), falls nötig.</li> <li>Stellen Sie sicher, dass die Messzelle ausreichend äquilibriert ist. Stellen Sie sicher, dass Sie entgaste Lösungsmittel verwenden.</li> <li>Wird der Vorgang wiederholt abgebrochen, wenden Sie sich an den Kundendienst.</li> </ul>

Meldung und Code	Beschreibung und Abhilfe
Code 8111 - Code 8115 Calibration of xx failed.	Die Kalibrierung von xx ist fehlgeschlagen. Wobei x = Name der Komponente, für welche die Kalibrierung
	fehlgeschlagen ist
	Mögliche Ursachen
	Die Raman-Wellenlängenkalibrierung ist fehlgeschlagen.
	Abhilfen
	<ul> <li>Beachten Sie die Bedingungen für die Wellenlängenkalibrierung (siehe Durchführung einer Wellenlängen-Kalibrierung (▶ Seite 137)) und Wellenlängenvalidierung (siehe Durchführung einer Wellenlängenvalidierung (▶ Seite 139)).</li> </ul>
	<ul> <li>Prüfen Sie, dass die Messzelle korrekt eingebaut ist und dass die Schrauben fingerfest angezogen sind.</li> </ul>
	<ul> <li>Reinigen Sie die Messzelle (siehe Reinigen der Messzelle (&gt; Seite 144)), falls nötig.</li> </ul>
	<ul> <li>Stellen Sie sicher, dass die Messzelle ausreichend äquilibriert ist und keine Luftblasen vorhanden sind. Stellen Sie sicher, dass Sie entgaste Lösungsmittel verwenden.</li> </ul>
	Wird der Vorgang wiederholt abgebrochen, wenden Sie sich an den Kundendienst.
Code 8116 and 8118 to 8121	Fehler beim Initialisieren
Initialization failed	Mögliche Ursachen
	<ul> <li>Der Detektor konnte nicht initialisiert werden.</li> </ul>
	Eventuell ist die Messzelle kontaminiert.
	Abhilfen
	<ul> <li>Prüfen Sie, dass eine Messzelle korrekt eingebaut ist.</li> </ul>
	<ul> <li>Pr üfen Sie, dass die sich die Eluentenzusammensetzung in der Messzelle nicht  ändert und es keine Luftblasen gibt. Stellen Sie sicher, dass Sie entgaste Lösungsmittel verwenden.</li> </ul>
	<ul> <li>Reinigen Sie die Messzelle (siehe Reinigen der Messzelle (&gt; Seite 144)), falls nötig.</li> </ul>

# 8.3 Störungen beim Betrieb

Dieser Abschnitt gibt Ihnen einen Überblick über mögliche Störungen beim Betrieb und Abhilfen.

### 8.3.1 Beheben von Undichtigkeiten

Wann

Der Leaksensor ist feucht. Der Leaksensor hat eine Undichtigkeit erkannt.

Erforderliche Teile und zusätzliche Utensilien

- Ersatzteil, je nach Erfordernis
- Tuch oder Papiertuch

#### Vorbereitungen

Wenn Sie Undichtigkeiten beseitigen, beachten Sie die Sicherheitshinweise und allgemeinen Regeln für Wartung und Service in Wartung und Service (▶ Seite 127).

Gehen Sie wie folgt vor

1. Finden Sie die Ursache der Undichtigkeit. Undichtigkeiten treten in der Regel an Verbindungen auf. Undichtigkeiten können jedoch auch in einer Messzelle auftreten.

Position der Undichtigkeit	Schritte
Messzelleneinlass und/oder -auslass	1. Ziehen Sie die Verbindung nach, an der Flüssigkeit sichtbar ist.
	<ol> <li>Wenn die Verbindung dicht zu sein scheint, jedoch immer noch undicht ist, entfernen Sie die Kapillare/Fitting und prüfen Sie, diese auf Beschädigung.</li> <li>Ersetzen Sie die Einlasskapillare oder Waste-Leitung falls erforderlich.</li> </ol>
Leckage-Wanne, aber nicht Messzelleneinlass oder -auslass	<ol> <li>Entfernen Sie die Messzelle aus dem Detektor und pr</li></ol>
	2. Wenn an der Messzelle Anzeichen von Undichtigkeit zu erkennen sind, trocknen Sie den Messzellenschacht vorsichtig und lassen Sie Restfeuchte verdunsten, bevor Sie die Messzelle ersetzen.

- 3. Saugen Sie mit einem Tuch oder Papiertuch alle Flüssigkeit gründlich auf, die sich in der Leckage-Wanne und unter dem Leaksensor angesammelt hat. Achten Sie darauf, den Sensor nicht zu verbiegen.
- 4. Lassen Sie dem Sensor einige Minuten Zeit, sich auf die Umgebungstemperatur einzustellen.
- 5. Wenn der Leaksensor keine Undichtigkeit mehr meldet, können Sie den Betrieb wieder aufnehmen.

#### Sehen Sie dazu auch

- Ausbauen der Messzelle ( Seite 142)
- Einbauen der Messzelle (
  Seite 145)

#### 8.3.2 Weitere mögliche Gerätestörungen beim Betrieb

In diesem Abschnitt finden Sie Informationen zu weiteren Störungen, die beim Betrieb des Vanquish-Gerätes auftreten können. Suchen Sie in der Tabelle, welche die entsprechende Art von Fehlerbild behandelt, nach möglichen Ursachen und lesen Sie die Beschreibung zur Abhilfe, um das Problem schnell zu lösen.

Prüfen Sie auch den Chromeleon Audit Trail auf eine entsprechende Meldung, wenn eine Störung auftritt. Die Meldung enthält gegebenenfalls weitere Informationen.

Beachten Sie, dass dieser Abschnitt nur Informationen zu Fehlerbildern und Ursachen liefert, die direkt mit dem Vanquish-Gerät zusammenhängen. Weitere Informationen zur Fehlersuche im Vanquish-System finden Sie in der Vanquish-System-Betriebsanleitung.

**TIPP** Wenn Sie eine Störung mithilfe der Beschreibungen in diesem Kapitel nicht selbst beheben können, oder Sie auf Probleme stoßen, die hier nicht beschrieben sind, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst, um Unterstützung zu erhalten.

Störung	Mögliche Ursache	Abhilfe
Peak-Tailing	Einlass und Auslass der Messzelle in einem Fluoreszenz-Detektor ausgetauscht	Prüfen Sie, dass der Einlass und Auslass der Messzelle korrekt angeschlossen sind.
Negative Peaks	Die Absorption/ Fluoreszenz des Analyts ist geringer als die der mobilen Phase.	Wählen Sie eine andere UV- oder Fluoreszenz- Detektionswellenlänge aus. Verwenden Sie eine mobile Phase mit geringerer Absorption/Fluoreszenz. Lösen Sie die Probe in der mobilen Phase.

Störung	Mögliche Ursache	Abhilfe
	Die Fluoreszenz der Substanz oder des Eluents wird durch andere Bestandteile gequencht.	Gegebenenfalls können die negativen Peaks zur Auswertung herangezogen werden.
Linien im Fluoreszenzsignal	Die Empfindlichkeit wurde automatisch reduziert	Wählen Sie eine niedrigere Empfindlichkeit für Ihre Anwendung. Der Audit Trail informiert Sie, zu welcher Empfindlichkeit reduziert wurde.
Spikes	Xenon-Blitzlampe veraltet, defekt, oder nicht richtig installiert	Wenden Sie sich an den Kundendienst.
	Die Messzellentemper- atur des Fluoreszenz- Detektors liegt nahe dem Siedepunkt der mobilen Phase.	Schalten Sie die Temperaturregelung für die Messzelle aus oder verringern Sie die Einstellung der Temperatur.
Ungeeignete Geräte- einstellungen	Ungeeignete Wellenlänge, z.B. auf einer Flanke des UV- Spektrums	Wählen Sie eine Wellenlänge oder ein Anregung/ Emissions-Wellenlängenpaar nahe des Spektrenmaximums.
	Die Empfindlichkeit des Fluoreszenz-Detektors wurde automatisch reduziert.	Wählen Sie eine niedrigere Empfindlichkeit für Ihre Anwendung.
Starke Basisliniendrift	Die Absorption des Eluenten verändert sich, wenn der Gradient gefahren wird	Absorbierende Additive können das Absorptionsspektrum verändern, abhängig vom Lösungsmittel. Erwägen Sie variierende Additivkonzentrationen, um die Drift auszugleichen.
Unregelmäßige Schwankungen der Basislinie, starkes Rauschen	Xenon-Blitzlampe im Fluoreszenz-Detektor zu alt	Wenden Sie sich an den Kundendienst.
Messzelle im Fluoreszenz- Detektor erreicht nicht die eingegebene Temperatur	Messzellentemperatur zu hoch oder zu niedrig eingestellt	Bei hohen Flüssen kann die Messzelle gegebenenfalls die gewünschte Temperatur nicht erreichen. Wählen Sie eine niedrigere Messzellentemperatur.
		Die Temperatur kann aber auch zu niedrig gewählt worden sein, zum Beispiel unter der Raumtemperatur. Passen Sie die Einstellung der Temperatur an oder schalten Sie die Temperaturregelung aus.

# 9 Spezifikationen

In diesem Kapitel finden Sie die technische Spezifikation und die Leistungsspezifikation sowie Informationen zu den Materialien, die im Flussweg des Gerätes verwendet werden.

# 9.1 Leistungsspezifikationen

### 9.1.1 Detektor

### Die Leistungsdaten des Gerätes sind wie folgt spezifiziert:

Art	Spezifikation
Optisches Design	Zwei Monochromatoren mit holographischen, konkaven Gittern und elliptischen Spiegeln für eine hoch effiziente Lichtübertragung
Lichtquelle	Xenon-Blitzlampe (15W); Es kann zwischen drei verschiedenen Blitzfrequenzen (HighPower, Standard, LongLife) gewählt werden.
Referenzsystem	Referenzsensor hinter der Messzelle zum Ausgleich von Fluktuationen der Lampenintensität
Wellenlängenbereich	VC-D50: Anregung: 200 nm - 630 nm Emission: 265 nm - 650 nm Anzahl der Photomultiplier: 1 VC-D51: Anregung: 200 nm - 880 nm Emission: 265 nm - 900 nm Anzahl der Photomultiplier: 2 VF-D50: Anregung: 200 nm - 630 nm Emission: 220 nm - 650 nm Anzahl der Photomultiplier: 1 VF-D51: Anregung: 200 nm - 880 nm Emission: 220 nm - 900 nm Anzahl der Photomultiplier: 2
Spektrale Bandbreite	Anregung: 20 nm Emission: 20 nm
Spektrenscanmodi	Single Spectrum Scan Modi: Excitation, Emission, Synchronous Unter der Software Chromeleon 7: FL Field Acquisition Scan Modi: Excitation, Emission, Synchro
Emissionsfilter	VC-D50, VC-D51: fester Filter mit 280 nm VF-D50, VF-D51: Fünf programmierbare Positionen (Offen, 280 nm, 370 nm, 435 nm, 530 nm)
Schaltzeit Anregungs-/ Emissionswellenlänge	< 250 ms
Anzahl der Kanäle	VC-D50, VC-D51: ein Kanal VF-D50, VF-D51: bis zu 4 Kanäle

Art	Spezifikation
Datenaufnahmerate	Einstellbar: Bis zu 100 Hz (VC-D50, VC-D51) Bis zu 200 Hz (VF-D50, VF-D51 unter der Software Chromeleon 7 im Single- Channel-Modus (Einkanal-Modus)) Bis zu 4 Hz (VF-D50, VF-D51 im Multi-Channel-Modus (Mehrkanal-Modus))
Empfindlichkeit	Bis zu 4 Hz (VF-D50, VF-D51 im Multi-Channel-Modus (Mehrkanal-Modus)) PMT 1: Raman Signal-Rausch-Verhältnis: >550 ASTM über die gesamte Lebensdauer der Lampe; Testbedingungen: Standard-Messzelle, Wasser in Fluoreszenzqualität, Anregung 350 nm, Emission 397 nm (>2100 mit Dunkelstrom als Referenz); Testbedingungen: Standard- Messzelle, Wasser in Fluoreszenzqualität, Anregung 350 nm, Emission 450 nm PMT 2 (nur VF-D51): Raman Signal-Rausch-Verhältnis: >225 ASTM; Testbedingungen: Standard-Messzelle, Wasser in Eluoreszenzqualität, Anregung 350 nm, Emission 397 nm
	<ul> <li>(&gt;1050 mit Dunkelstrom als Referenz); Testbedingungen: Standard-</li> <li>Messzelle, Wasser in Fluoreszenzqualität, Anregung 350 nm, Emission</li> <li>450 nm</li> </ul>
Wellenlängengenauigkeit	± 2 nm (über die Detektor-Lebensdauer; einzeln für Anregungs- und Emissions-Monochromatoren)
Wellenlängenwiederholbarkeit	± 0,2 nm
Wellenlängenkalibrierung	Interne Kalibrierung, Anregungs-Monochromator mit Emissionslinien der Xenon-Blitzlampe, Emissions-Monochromator mit Raman-Verschiebung von Wasser und Emissionslinien der Xenon-Blitzlampe.
Wellenlängenvalidierung	<ul> <li>Interne Validierung, Anregungs-Monochromator mit Emissionslinien der Xenon-Blitzlampe, Emissions-Monochromator mit Raman-Verschiebung von Wasser und Emissionslinien der Xenon-Blitzlampe.</li> <li>Da die Wellenlängenbestimmung der Ramanbande sowohl von der Anregungs- als auch von der Emissionswellenlängengenauigkeit abhängig ist, ist die erlaubte Abweichung ± 3 nm.</li> </ul>
USB	1 USB-Port (USB 2.0, Typ "B") 1 USB-Hub mit 3 Ports (USB 2.0, Typ "A")
E/A-Schnittstelle	2 digitale E/A-Ports (Mini-DIN), jeweils mit einem digitalen Eingang und einem Relaisausgang
System Interlink	2 System Interlink -Ports (RJ45-8)
Analogausgang	2 Analogausgänge über optionale Erweiterungskarte zur Ausgabe von Emissionskanälen Auflösung: 20 bit Maximale Datenrate: 50 Hz Ausgänge über Software einstellbar (Ausgangsspannungsbereich 0 bis 1 V oder 0 bis 10 V, Empfindlichkeit und Offset)

Art	Spezifikation
Steuerung	Chromeleon 7
	Das Gerät kann auch mit anderen Datensystemen betrieben werden. Weitere Fragen beantwortet Ihnen gern die Thermo Fisher Scientific- Vertriebsorganisation.
	Tastatur mit 2 Funktionstasten zur Ausführung bestimmter Funktionen direkt am Gerät
Materialien im Flussweg	Siehe Spezifikationen für die Messzellen.
Sicherheitsmerkmale	Überprüfung der Optik und der Motoren beim Einschalten (Initialisierung) Kontrolle der Kühlventilatoren und der Elektronik Leakerkennung und sichere Leakbehandlung Identifizierung und Dokumentation des Messzellentyps
Good Laboratory Practice (GLP)	Alle Systemparameter werden im Chromeleon Audit Trail protokolliert. Funktionen zur Kontrolle der Betriebs- und Nutzungsbedingungen des Geräts. Dies beinhaltet die Überprüfung des Lampenalters, die Arbeitslast des PMT, des Gitters und der Filterbewegungen, sowie Service- und Qualifikationsintervalle über die Software Chromeleon.

## 9.1.2 Messzellen

Art	Standard-Messzelle	Mikro-Messzelle	
Messzellenvolumen	8 μL	2 μL	
Volumen Wärmetauscher und/oder Einlasskapillare	6,3 μL	3,3 μL	
Druckgrenze	2 MPa	4 MPa	
Verwendungs- empfehlungen	Größere Empfindlichkeit und ein besseres Signal-Rausch-Verhältnis, speziell für Säulen mit > 2,1 mm ID	Beste Auflösung in der UHPLC, speziell für Säulen mit ≤ 2,1 mm ID	
Temperatur- regelung	15 °C über Raumtemperatur bis 50 °C absolut		
Biokompatibilität	ja		
Materialien im Flussweg	Quarzglas, MP35N, Peek, Titan, Fluorpolymere ACHTUNG Informationen zur chemischen Beständigkeit der Materialien können Sie der technischen Literatur entnehmen.		
Informationen zu Lösungsmitteln und Additiven	Siehe Informationen zu Lösungsmitteln und Additiven (▶ Seite 27). Maximal erlaubte Lösungsmitteltemperatur: 80 °C.		
Good Laboratory Practice (GLP)	Identifikationschip		

### Die Leistung der Messzellen ist wie folgt spezifiziert:

# 9.2 Technische Spezifikationen

Die technischen Daten des Gerätes sind wie folgt spezifiziert:

Art	Spezifikation
Verwendungsbereich	Ausschließlich im Innenbereich
Umgebungstemperatur (Betrieb)	5 °C - 35 °C
Umgebungstemperatur (Lagerung)	-20 °C - 45 °C
Umgebungsluftfeuchtigkeit (Betrieb)	20% - 80% relative Feuchte (nicht-kondensierend)
Umgebungsluftfeuchtigkeit (Lagerung)	Maximal 60% relative Feuchte (nicht-kondensierend)
Betriebshöhe	Maximal 2000 m über Normalnull
Verschmutzungsgrad	2
Leistungsaufnahme	Weiter Bereich, 100 – 240 V AC, ± 10 %; 50/60 Hz; max. 245 W / 255 VA
Überspannungskategorie	Ш
Emissionsschalldruckpegel	< 50 dB (A)
Abmessungen (Höhe x Breite x Tiefe)	15,9 x 42 x 62 cm
Gewicht	Circa 21 kg
# 10 Zubehör, Ersatzteile und Verbrauchsmaterialien

In diesem Kapitel ist das Standard-Zubehör aufgeführt, das zusammen mit dem Gerät ausgeliefert wird, sowie Zubehör, das optional bestellt werden kann. Zusätzlich finden Sie Informationen zur Nachbestellung von Verbrauchsmaterialien und Ersatzteilen.

# **10.1** Allgemeine Informationen

Das Gerät darf ausschließlich mit Ersatzteilen und zusätzlichen Komponenten, Optionen und Peripheriegeräten betrieben werden, die von Thermo Fisher Scientific ausdrücklich autorisiert und freigegeben sind.

Zubehör, Ersatzteile und Verbrauchsmaterialien werden laufend dem neuesten technischen Stand angepasst. Eine Änderung der Bestellnummern ist deshalb nicht auszuschließen. Wenn nicht anders angegeben, werden jedoch bei Bestellung der aufgeführten Bestellnummern stets voll kompatible Teile geliefert.

# 10.2 Zubehörkit

Das Zubehörkit enthält die in der Tabelle aufgeführten Teile. Der Inhalt des Zubehörs kann jederzeit geändert werden und von den in dieser Anleitung enthaltenen Angaben abweichen. Aktuelle Informationen zum Inhalt des Zubehörkits bei Auslieferung des Geräts finden Sie in der beiliegenden Zubehörliste.

Zubehörkit

Artikel	Menge im Zubehör
Einsatz für die Zwischenwand, um isolierte Kapillaren zu führen	1
System-Interlink-Kabel	1
Schlauchklammer	1
USB-Kabel, USB 2.0, High-speed, Typ A auf Typ B	1
Viper-Kapillare, Säulenthermostat - Fluoreszenz-Detektor	1
Viper-Kapillare, UV-/VIS-Detektor - Fluoreszenz-Detektor, 0.18 mm I.D.	1
Waste-Leitung	1

Informationen zur Nachbestellung finden Sie unter Ersatzteile und Verbrauchsmaterialien (▶ Seite 185).

# 10.3 Optionales Zubehör

Messzellen und Messzellenzubehör

Artikel	BestNr.
Standard-Messzelle, biokompatibel	6079.4230
Mikro-Messzelle, biokompatibel	6079.4330
Spül- und Injektionskit mit Spritze für Messzellen, einschließlich Spritze	6078.4200
Viper-Kapillare, I.D. x Länge 0,18 x 450 mm, Edelstahl Zur Verbindung der Säule mit dem Messzelleneinlass, wenn eine UltiMate 3000 Messzelle verwendet wird.	6040.2365
Überdruck-Entlüftungsventil Geeignet zum Schutz der Mikro-Messzelle im Fluoreszenz- Detektor gegen Überdruck. Das Ventil öffnet bei einem Druck von 4 MPa (40 bar).	6079.9240

**TIPP** Messzellen von UltiMate 3000-Fluoreszenz-Detektoren haben den Einlass und Auslass auf der linken Seite. Um den Messzellenauslass an der linken Seite der Messzelle in einem Vanquish-Fluoreszenz-Detektor anzuschließen, empfehlen wir die Verwendung der 450 mm langen Kapillare, die in der Tabelle weiter oben genannt ist.

#### Sonstiges

Artikel	BestNr.
DAC-Einschub Stellt zwei Analogausgänge zur Verfügung. Wenden Sie sich für den Einbau an den technischen Kundendienst von Thermo Fisher Scientific.	6083.0900
Dual-PMT-Option Enthält einen zweiten PMT für den nahen Infrarotbereich zum Nachrüsten. Wenden Sie sich für den Einbau an den technischen Kundendienst von Thermo Fisher Scientific.	6078.5360

# 10.4 Ersatzteile und Verbrauchsmaterialien

#### Schläuche und Kapillaren

Beschreibung	BestNr.	
Viper-Kapillare, 350 mm Länge, MP35N, zur Verbindung zum Säulenthermostaten	6042.2340	
Viper-Kapillare, 300mm Länge, MP35N, zur Verbindung von einem UV-/VIS-Detektor zum Fluoreszenz-Detektor	6042.2322	
Waste-Leitung	6036.2425	
Verschlussstopfen für den Einlass und Auslass der Messzelle	6200.5502	
Systemkapillaren, siehe Vanquish-System-Betriebsanleitung.		

#### Sonstiges

Beschreibung	BestNr.
Gerätetüren-Kit mit rechter Tür und linker Tür	6083.3018
Sicherungskit, Vanquish-System Das Kit enthält die passenden Sicherungen für die Vanquish- Systemmodule. Verwenden Sie für den Detektor ausschließlich Sicherungen des Typs 3.15 AT, 250 V AC, träge.	6036.0002
Verpackungsmaterial für Detektor mit Gehäuse	6083.0090

#### Schnittstellen-Kabel

Beschreibung	BestNr.
Digitales I/O-Signalkabel, 6-polig, Kabellänge: 5 m	6036.0006
System-Interlink-Kabel	6036.0004
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0 Kabellänge: 0,5 m	6720.8910
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0 Kabellänge: 1 m	6035.9035
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0 Kabellänge: 5 m	6911.0002

#### Netzkabel

Beschreibung	BestNr.
Netzkabel, Australien	6000.1060
Netzkabel, China	6000.1080
Netzkabel, Dänemark	6000.1070
Netzkabel, EU	6000.1000
Netzkabel, Indien, SA	6000.1090

Beschreibung	BestNr.
Netzkabel, Italien	6000.1040
Netzkabel, Japan	6000.1050
Netzkabel, UK	6000.1020
Netzkabel, US	6000.1001
Netzkabel, Schweiz	6000.1030

# 11 Anhang

In diesem Kapitel finden Sie weitere Informationen zur Konformität, zu den UV-Cutoff-Wellenlängen und zur Verwendung der digitalen E/A-Schnittstellen.

# 11.1 Informationen zur Konformität

#### 11.1.1 Konformitätserklärungen

#### CE-Konformitätserklärung

Das Gerät entspricht den Anforderungen für die CE-Kennzeichnung und genügt den geltenden Anforderungen.

#### cTUVus-Konformität

Das cTUVus-Zeichen auf dem Gerät zeigt an, dass das Gerät die Anforderungen für das cTUVus-Zeichen erfüllt. Die Erfüllung der geltenden Anforderungen wurde durch die TÜV Rheinland of North America Inc. geprüft.

#### RoHS-Konformität

Dieses Produkt entspricht den Richtlinien für RoHS (Restrictions of Hazardous Substances):

 Europäische RoHS-Richtlinie
 Richtlinie zur Beschränkung der Verwendung bestimmter gefährlicher Stoffe in elektrischen und elektronischen Geräten

Das CE-Zeichen auf dem Gerät gibt an, dass das Gerät die Vorgaben der Richtlinie erfüllt.

 China-RoHS-Richtlinien Measures for Administration of the Pollution Control of Electronic Information Products (Maßnahmen zur Kontrolle von Umweltverschmutzungen durch elektronische Produkte)

Folgende Logos können sich auf dem Gerät befinden:

Logo	Beschreibung
Das grüne Logo kennzeichnet Geräte, die keine in den Richtl genannten gefährlichen Stoffe enthalten.	
	Das orangene Logo mit einer ein- oder zweistelligen Zahl kennzeichnet Geräte, die in den Richtlinien genannte gefährliche Stoffe enthalten. Die Zahl gibt den EFUP-Zeitraum (Environment- Friendly Use Period, Zeitraum, in dem die umweltfreundliche Nutzung gegeben ist) an. Das Gerät verursacht (bei bestimmungsgemäßer Verwendung) während dieses Zeitraums keine Schäden für die menschliche Gesundheit oder die Umwelt.
	Weitere Informationen finden Sie auf http://www.thermofisher.com/ us/en/home/technical-resources/rohs-certificates.html

#### 11.1.2 WEEE-Konformität

Dieses Produkt erfüllt die Bestimmungen der EU-Richtlinie über Elektround Elektronik-Altgeräte (WEEE-Richtlinie). Es ist mit dem folgenden Symbol gekennzeichnet:



#### Abbildung 43: WEEE-Symbol

Thermo Fisher Scientific hat in jedem Europäischen Unions-Mitgliedstaat (EU-Mitgliedstaat) Verträge mit einem oder mehreren Wiederverwertungs- oder Entsorgungsunternehmen abgeschlossen; dieses Produkt sollte zur Entsorgung oder Wiederverwendung an diese Partner übergeben werden. Weitere Fragen beantwortet Ihnen Thermo Fisher Scientific gern.

#### 11.1.3 Einhaltung der FCC-Richtlinien

Dieses Gerät wurde geprüft und erfüllt die Grenzwerte für Digitalgeräte der Klasse A gemäß Absatz 15 der amerikanischen FCC-Richtlinien.

Diese Grenzwerte sollen einen angemessenen Schutz gegen elektromagnetische Störungen beim Betrieb in gewerblich genutzten Räumen gewährleisten. Das Gerät erzeugt und verwendet Hochfrequenzenergie und kann diese auch selbst aussenden. Bei nicht ordnungsgemäßer Installation und Verwendung gemäß der Betriebsanleitung sind schädliche Störungen des Funkverkehrs möglich.

#### 11.1.4 Versionsgeschichte der Anleitung

Version	In der Anleitung beschrieben	
2.0	VC-D50, VC-D51, VF-D50, VF-D51	
1.2a	VF-D50, VF-D51	
1.2	VF-D50, VF-D51	
1.1	VF-D50, VF-D51	

Die Anleitung wurde auf Englisch erstellt (Originalanleitung). Andere Sprachversionen sind Übersetzungen der englischen Originalanleitung.

# 11.2 UV-Cutoff-Wellenlängen von Lösungsmitteln

Die UV-Cutoff-Wellenlänge ist die niedrigste verwendbare Messwellenlänge. Die Zusammensetzung der mobilen Phase wirkt sich auf die UV-Cutoff-Wellenlänge aus.

Im Allgemeinen bestehen mobile Phasen aus Lösungsmitteln wie Wasser, Acetonitril, Methanol oder anderen Stoffen. Es können auch Salze, wie Natriumhydroxid (NaOH), enthalten sein.

Die meisten Lösungsmittel haben eine UV-Cutoff-Wellenlänge innerhalb des spektralen Bereichs des Geräts. Führen Sie für optimale Messergebnisse quantitative Messungen bei einer Wellenlänge durch, die ausreichend über der UV-Cutoff-Wellenlänge liegt. Die UV-Cutoff-Wellenlänge hängt auch von der Qualität des Lösungsmittels ab. Um Informationen zur UV-Cutoff-Wellenlänge für ein Lösungsmittel zu erhalten, wenden Sie sich an den Hersteller des Lösungsmittels.

Der Brechungsindex von verschiedenen Lösungsmitteln kann unterschiedlich sein. Das Verändern der Zusammensetzung der mobilen Phase kann daher zu Basislinienschwankungen führen.

# 11.3 Digital I/O

Die digitalen I/O-Ports (Dig I/O) können zur Übertragung von digitalen Signalen an externe Geräte verwendet werden. Jeder Port verfügt über:

- einen digitalen Eingang
- einen Relaisausgang

Anschlussbelegung



Abbildung 44: Digital I/O-Port

Pin	Beschreibung
1	Nicht verwendet
2	Relaisausgang — Relay_NC (NC = Normally Closed = Ruhekontakt)
3	Ground — GND
4	Digitaleingang — Input
5	Relaisausgang — Relay_COM COM ist der gemeinsame Kontakt für NO und NC. Ist das Relais nicht aktiviert oder ist das Gerät ausgeschaltet, besteht eine Verbindung zwischen COM und NC. Ist das Relais aktiviert, besteht eine Verbindung zwischen COM und NO.
6	Relaisausgang — Relay_NO (NO = Normally Open = Arbeitskontakt)

In der folgenden Tabelle finden Sie die Zuordnung der Funktionen zu den Steckerpins und zur Farbe der Kabelader, die mit dem jeweiligen Pin verbunden ist.

Pin	Farbe Ader	Signalname	Signalpegel	Bemerkung
1	Rosa			Nicht verwendet
2	Grau	Relaisausgang — Relay_NC	Potentialfrei 0-24 V, 0-100 mA	Öffner
3	Grün	Ground — GND	Ground	Bezugspotential
4	Gelb	Digital input — Input	Eingang (low active): Ein: 0-0.4 V Aus: 2.2-5 V Pullup-Widerstand: 47 kΩ bis 5 V	<ul> <li>Digitaleingang; das</li> <li>Referenzpotential ist Ground.</li> <li>Beachten Sie folgende Hinweise:</li> <li>Die maximale Eingangsspannung darf am Eingang +5 V in Bezug auf Ground nicht überschreiten.</li> <li>Die minimale Eingangsspannung darf das Groundpotential nicht unterschreiten.</li> </ul>
5	Weiß	Relaisausgang — Relay_COM	Potentialfrei	Mittelkontakt für NO und NC
6	Braun	Relaisausgang — Relay_NO	Potentialfrei 0-24 V, 0-100 mA	Schließer

#### Voraussetzungen

Um die Digital-I/O-Funktionalität nutzen zu können, müssen folgende Voraussetzungen erfüllt sein:

- Der digitale I/O-Port muss mit dem externen Gerät über das digitale I/O-Signalkabel (Best.-Nr. 6036.0006) verbunden sein.
- Die Eingänge und Ausgänge, die Sie verwenden möchten, müssen im Instrument Configuration Manager ausgewählt sein.

#### Verbinden des Digitalen I/O-Ports

- 1. Stecken Sie den 6-poligen Stecker des Kabels in den digitalen I/O-Port, den Sie verwenden möchten.
- 2. Verbinden Sie für jeden Relaisausgang oder digitalen Eingang, den Sie verwenden möchten, die geeignete Signalader und Groundader mit dem entsprechenden Anschluss am externen Gerät. Einzelheiten entnehmen Sie bitte der Dokumentation für das externe Gerät.

Auswählen der Eingänge und Ausgänge in der Software Chromeleon

 Wählen Sie im Dialog für das Gerät, auf den Seiten Inputs und Outputs, die Eingänge und Ausgänge aus, die Sie verwenden möchten. Die Nummerierung im Dialog entspricht der Nummerierung am Port.

# Index

#### Α

#### B

Basislinienrauschen	109
Basislinienspektrum	104
Basislinienverhalten	109, 120
Bestellinformationen	181
Betrieb	41, 91
Bedienelemente	
Messzellen	100
Sicherheitshinweise	
Stromversorgung ein-/ausschalten	
Unterbrechen	121
wichtige Parameter	106
Betriebsmodi	101

## С

CE-Kennzeichnung	188
Chloridkonzentration	28
Chromeleon	41
Audit Trail	164
Einstellen der Wellenlängen	110
Geräteeinrichtung	89
mehrere Kanäle messen	102
Moduleinrichtung	89

106
135
121
121
99
188

## D

Datenaufnahmerate 1	109, 117
Hinweise	117
Dekontaminieren	133
Dekontaminierung	133
Detektor	
Empfindlichkeit	111
Weitere mögliche Störungen beim Bet	rieb
	172
Detektor-Waste-Leitung	76
Dig I/O	58, 191
Digital I/O	58, 191
Drainage	77

## Ε

Einlasskapillare		72
Einrichtung		
Hardware		54
Software		89
Systemaufbau		54
Einschalten		88
Einstellungen	1	106
Emission Scan	1	104
Emissionswellenlänge	109, 1	L10
Empfindlichkeit	109, 1	111
erlaubter Emissionswellenlängenbereich	1	109
Ersatzteile	181, 1	185
Excitation Scan	1	104

#### F

FCC	189
Fehlersuche	163
Allgemeine Informationen	164
Meldungen	166
Störungen beim Betrieb	171
Filterrad 109,	115
Firmware-Update	148

FL Field Acquisition	. 105
Fluoreszenzintensität	. 109
Führungsloch	65
Funktionsprinzip	33

## G

Gerät	
Außerbetriebnahme	121
Neustart nach langfristiger	
Außerbetriebnahme	125
Optimierung	108
Vorbereiten für Betrieb	. 98
Gerätestart nach langfristiger	
Außerbetriebnahme	125
Gesetzliche Bestimmungen	. 29

#### Н

Handschuhe	23
HighPower-Modus	119
Hinweise	
Betrieb	93
Installation	48
Messzellen Handhabung	100
Wartung	129
Wartung der Messzelle	141

#### I

Identifikationschip	37
Messzelle	37
Inbetriebnahme	98
Initialize	107
INIT-Taste	95
Innenansicht	36
Installation	47
Anforderungen an den Aufstellungsort	51
Kapillaren und Schläuche	65
Sicherheitshinweise	48
System	49
Interlink	57

#### Κ

Kalibrierung 1	37
Kapillaren	
Führung	65
Installieren	69
Viper	69

Kondensation	52,	61
Kurzzeitige Außerbetriebnahme	1	21

#### L

Lampe	39 <i>,</i> 136
Lampen-Modus	119
Tauschen	136
Lampenalter	136
Lampenlebensdauer	136
Lampen-Modus	109
Langfristige Außerbetriebnahme	122
Langpassfilter	115
Leakerkennung	40, 107
Leaksensor	171
Leaksensor-Einstellung	107
LED-Leiste	94, 95 <i>,</i> 164
Lieferumfang	46
LongLife-Modus	120
Lösungsmittel	
Chloridkonzentration	28
Information	27
pH-Bereich	27

#### Μ

MaxPMTSaturation	112
Mehrfach-Wellenlängenbetrieb	101, 102
Meldungen	166
Messzelle	37
Anschluss des Auslasses	76
Anschluss des Einlasses	72
Entfernen	142
Erstinstallation	62
Flussverbindungen	71
Handhabung	141
Hinweise	100
Identifikationschip	37
Installation	145
Reinigen	144
Rückdruckermittlung	78
Tauschen	142, 145
Temperaturregelung	38
Typen	38
Messzellen-Temperatur	109, 118
Methodenentwicklung	108
Moduleinschub	156
Entfernen	156
Montieren	159
Zurückschicken	158

Moduleinschub entfernen	156
Moduleinschub installieren	159
Multi-Channel Performance	102
Multi-Channel-Modus	101
Mute Alarm	. 94

#### Ν

Nachweisgrenze	109
Netzkabel	52
Nicht initialisiert	137

## 0

optimale Wellenlänge	110
Optimierung	108
Datenaufnahmerate	117
Hinweise	108
Parameter-Übersicht	109

### Ρ

Peakbreite	109
pH-Bereich	27
PMT	109, 116
Predictive Performance	135
Puffer	27
Information	27
Konzentration	27

## R

Reinigen	133
Reproduzierbarkeit	109
RoHS-Kennzeichnung	188

## S

Scans	
Einzelspektrum	103
FL Field Acquisition	105
Schläuche	69
Schlauchführungen	65
Schlauchkanal	65
Schlauchklammer	65
Schutzbrille	23
Schutzklasse	21
Schutzkleidung	23
Selektivität	109
Service	127

Sicherheitshinweise	
allgemein	21
allgemeine Risiken	24
Betrieb	
elektrische Sicherheit	23
Installation	48
Netzkabel	52
Qualifikation des Personals	22
Schutzausrüstung	22
Service	129
Verhalten im Notfall	
Wartung	129
Sicherheitssymbole	18, 19
Sicherungen	146
Signalwörter	18
Single Spectrum Scan (Erfassung von	
Einzelspektren)	103
Single-Channel-Modus	101
Smart Shutdown	121
Smart Standby	121
Smart Startup	
Speicherplatz	109
Spektrenscans	103
Spezifikationen	175
Leistung (Detektor)	176
Leistung (Messzellen)	178
technisch	179
Standard-Modus	120
STATUS LED	164
Statusanzeige	
LED-Leiste	94, 95, 164
STATUS LED	164
Status-LED	94, 95
Status-LED	94, 95
Störungen beim Betrieb	171
Stromversorgung (Hinweise)	51, 52
Stromversorgung ein-/ausschalten	
Synchronous Scan	104
System-Äquilibrierung	
Systemaufbau	54
Systemdrainage	77
System-Interlink	57

## Т

Tastatur	
Temperaturregelung	38
Transport	152
Transportieren	50

Tür	53
Aushängen	128, 150
Entfernen	128
öffnen	53
Tauschen	150
Typenschild	19

## U

Überblick (Funktionen)	31
Undichtigkeit	171
Undichtigkeiten	171
Universal Serial Bus	58
USB	58
UV-Cutoff-Wellenlängen	190

#### V

Verbrauchsmaterialien	181, 185
Versand	152
Verwendungszweck	20
Viper-Fitting System	69
Vorbereitung	
Modul entfernen	156

#### W

Wartung 127,	132
Allgemeine Regeln	131
Dekontaminieren	133
Einführung	128
Firmware-Update	148
Intervall	132
Lampe	136
Messzellen	141
Predictive Performance	135
Reinigen	133
Sicherheitshinweise	129
Sicherungen	146
Türen	150
Wellenlängenkalibrierung	137
Wellenlängenvalidierung	139
Wellness	135
Waste-Leitung	. 76
WEEE	189
Weitere mögliche Störungen beim	
Detektorbetrieb	172
Wellenlängen bestimmen	110
Wellenlängenkalibrierung	137
Wellenlängenschaltung	110

Wellenlängenvalidierung ..... 139

# Ζ

	102
46,	181
	184
	183
46,	183
	116
	 46,  46,

www.thermofisher.com

Thermo Fisher Scientific Inc. 168 Third Avenue Waltham Massachusetts 02451 USA

