

thermoscientific

Vanquish

Neo System VN-A10, VN-C10, VN-P10, VN-S10

Betriebsanleitung

4822.5001-DE Version 2.0 •

September 2024





Copyright © 2024 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

Übersetzung der Original-Betriebsanleitung

Die Hardware-Beschreibungen in dieser Anleitung beziehen sich auf die Gerätetypen: VN-A10-A-02, VN-C10-A-01, VN-P10-A-01, VN-S10-A-01.

Warenzeichen

Alle Warenzeichen sind Eigentum von Thermo Fisher Scientific und ihren Tochtergesellschaften sofern nicht anders angegeben.

Acrobat, Adobe, und Adobe Reader sind Warenzeichen oder eingetragene Warenzeichen von Adobe Systems Incorporated.

Microsoft und Windows sind Warenzeichen von Microsoft Corporation.

MP35N ist ein Warenzeichen von SPS Technologies.

PharMed ist ein Warenzeichen von Saint-Gobain Performance Plastics.

Torx ist ein Warenzeichen von Acument Intellectual Properties, LLC.

VICI ist ein Warenzeichen von Valco Instruments Co. Inc. und VICI AG International.

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_2 -Emissionen.

Kontaktinformationen

Wenn Sie Unterstützung von Thermo Fisher Scientific benötigen, finden Sie Informationen dazu auf der folgenden Webseite: https://www.thermofisher.com/us/en/home/technical-resources/ contact-us.html

Inhalt

1	Ver	wend	ung dieser Anleitung	15
	1.1	Über (diese Anleitung	16
	1.2	Konve	entionen für Sicherheitshinweise	17
	1.3	Beson	dere Hinweise	18
	1.4	Турод	rafische Konventionen	19
	1.5	Refere	enzdokumentation	20
2	Sic	herhei	it .	21
	2 1	Sichor	hoitesumhala und Signalwärtar	
	2.1	2 1 1	Sicherheitssymbole und Signalwörter in dieser Anleitung	22
		2.1.1	Sicherheitssymbole und Signalworter in dieser Amertung	22
		2.1.2	Sicherheitssymbole am System	22
		2.1.5	Sichemeitssymbole am System	25
	2.2	2.1.4		25
	2.2	verwe	endungszweck	25
	2.3	Sicher	heitsmaßnahmen	26
		2.3.1	Allgemeine Sicherheitsinformationen	26
		2.3.2	Qualifikation des Personals	27
		2.3.3	Persönliche Schutzausrüstung	27
		2.3.4	Sicherheitsmaßnahmen beim Umgang mit elektrischen Geräten	28
		2.3.5	Allgemeine Restrisiken	29
		2.3.6	Verhalten im Notfall	31
	2.4	Inforn	nationen zur Konformität	32
3	Üb	ersich	t über das System	33
	3.1	Syster	nmerkmale	34
	3.2	Besch	reibung des Systems	37
	3.3	Svster	nsockel	38
		3.3.1	Komponenten des Systemsockels	38
		3.3.2	Systemsockelschloss	39
	3.4	Syster	n Controller	40
	3.5	Pump	e	41
		3.5.1	Innenansicht der Pumpe	41

	3.6	.6 Autosampler		44
		3.6.1	Innenansicht des Autosamplers	44
		3.6.2	Funktionsprinzip des Autosamplers	46
		3.6.3	Anschlussbelegung der Autosampler-Ventile	48
		3.6.4	Probenraum	49
	3.7	Säulen	thermostat	53
		3.7.1	Innenansicht des Säulenthermostaten	53
		3.7.2	Funktionsprinzip des Säulenthermostaten	54
		3.7.3	Übersicht der Säulenschaltventile	54
	3.8	Solven	t Rack	56
	3.9	Leaker	kennung	57
	3.10	Nadel-	Waschsystem (Übersicht)	58
	3.11	Hinter	spülung (Übersicht)	60
	3.12	Ein- ur	nd Ausschalten	62
	3.13	Betriel	o des Systems	64
	3.14	Purger	n (Übersicht)	65
				~ 7
4	Aus	раске	n	6/
	4.1	Beschä	idigte Verpackung, Mängel bei Ankunft des Geräts	68
	4.2	Lieferu	ımfang	69
	4.3	Auspa	cken des Systems	70
	4.4	Auspa	cken des Säulenthermostaten	72
	4.5	Entspe	erren des Systemsockels	73
5	Inst	allatio	n	75
	5 1	Sicharbaitshinwaisa für die Installation		
	5.1	Anform		70
	5.2			70 70
		522	Hinweise zur Stromversorgung	70 70
		5.2.3	Netzkabel	80
		5.2.4	Kondensation	81
		5.2.5	Betriebsbedingungen	81
	5.3	Aufste	llen des Systems	83
	5.4	Öffner	n und Schließen der Gerätetüren	85

	5.5	Installa	ation des Säulenthermostaten	86
		5.5.1	Einbau des Säulenthermostaten in den Systemturm	86
		5.5.2	Installieren des Vanquish-Schaltventils	88
		5.5.3	Installieren des VICI-Schaltventils	90
		5.5.4	Installieren des passiven Vorheizers	92
		5.5.5	Installieren von Säulen im Säulenthermostaten	94
	5.6	Entfer	nen der Transportsicherungen des Autosamplers	96
		5.6.1	Entfernen der Isolierabdeckung	98
		5.6.2	Einbauen der Isolierabdeckung	99
	5.7	Montie	eren des Displays	100
	5.8	Anschl	ießen der Kabel	107
		5.8.1	Übersicht der Anschlüsse an den Systemmodulen	107
		5.8.2	Übersicht der Anschlüsse am System Controller	108
		5.8.3	Anschließen der USB-Kabel	109
		5.8.4	Anschließen der Display-Kabel	111
		5.8.5	Anschließen der System-Interlink-Kabel	113
		5.8.6	Anschließen eines LAN-Kabels	115
		5.8.7	Anschließen der Netzkabel	115
		5.8.8	Verlegen der Signalkabel in den Kabelhalterungen	116
	5.9	Anschl	ießen eines Massenspektrometers	117
	5.10	Sperre	n des Systemsockels	118
	5.11	Anschl	ießen der Flussverbindungen	119
		5.11.1	Kapillar- und Schlauchführung durch das System	119
		5.11.2	Anschließen von Fittingen, Kapillaren und Schläuchen	120
		5.11.3	Anschließen der Waste-Leitungen	123
	5.12	Füllen	der Lösungsmittelbehälter	128
	5.13	Füllen	der Behälter für die Nadel-Waschflüssigkeit	130
	5.14	Füllen	des Behälters für die Hinterspülflüssigkeit	133
	5.15	Einsch	alten des Systems nach der Installation	135
	5.16	Integri	eren in ein Geräte-LAN	137
	5.17	Vorbei	reiten des Systems für den ersten Betrieb	138
6	Bet	rieb		41
	6.1	Sicherl	heitshinweise zum Betrieb	142
	6.2	Inform	ationen zu Lösungsmitteln und Additiven	144
		6.2.1	Allgemeine Kompatibilität	144
		6.2.2	Erlaubte pH-Bereiche	144

	6.2.3	Erlaubte Konzentrationen	145	
	6.2.4	Verwenden von Lösungsmitteln und Additiven	145	
6.3	Allgen	Allgemeine Hinweise zum Systembetrieb 148		
6.4	Einsch	alten des Systems	149	
6.5	Bedier	nelemente	150	
	6.5.1	Tastaturen	150	
	6.5.2	LED-Statusleisten	153	
6.6	Vanqu	ish User Interface	155	
	6.6.1	System Controller-Informationen	155	
	6.6.2	Systemstatus-Symbole	155	
	6.6.3	Navigationsleistensymbole und andere Symbole	156	
	6.6.4	Betriebszustand des Systems	156	
	6.6.5	Startfenster	156	
	6.6.6	Skriptfenster	157	
	6.6.7	Einstellungsfenster	157	
6.7	Anpas	sen des Displays	158	
	6.7.1	Mögliche Positionen des Displays	158	
	6.7.2	Verstellen des Displays von der vorderen Position in die seitliche Pos	ition 159	
6.8	Vorbe	reiten des Systems für den Betrieb	160	
	6.8.1	Spülen des Systems	160	
	6.8.2	Äquilibrieren des Systems	162	
	6.8.3	Beladen des Drehtellers	163	
	6.8.4	Empfehlungen zu Kappen und Verschlüssen von Probengefäßen	166	
6.9	Optim	ieren der Leistung	167	
	6.9.1	Abhängigkeit von verfügbarem Pumpendruck und Flussrate	167	
	6.9.2	Hinzufügen von kundenspezifischen Lösungsmitteln zur Kalibriertabe	elle	
		·····	168	
	6.9.3	Andern des Lösungsmitteltyps	170	
	6.9.4	Andern der Hinterspültlüssigkeit	170	
	6.9.5	Andern der Nadel-Waschflussigkeit	170	
	6.9.6	Durchfuhren des Nullpunktabgleichs für die Flusssensoren der Pump	e 1/1	
	6.9.7	Waschen der Probennadel	. 172	
	6.9.8	Thermostatisierung des Autosampier-Probenraums	. 1/3	
<i>.</i>	0.9.9		1/3	
6.10	Außer	betriebnenmen des Systems	175	
	6.10.1	Kurze Betriebsunterbrechung des Systems	175	
	6.10.2	Lange Betriebsunterbrechung des Systems	175	

7	Wa	rtung	und Service	L77
	7.1	Einfüh	rung in Wartung und Service	178
	7.2	Sicher	heitshinweise zu Wartung und Service	179
	7.3	Allgen	neine Regeln für Wartung und Service	182
	7.4	Wartu	ing und Wartungsintervalle	184
		7.4.1	Wartungszeitplan	184
		7.4.2	Spülen der Pumpe zur Vermeidung von Kontamination	187
		7.4.3	Spülen bei hartnäckiger Verschmutzung	189
		7.4.4	Reinigen oder Dekontaminieren eines Systemmoduls oder des System	S 100
	7.5	Wartu	ing des Pumpenkopfs	190
		7.5.1	Überblick über die Pumpenkopfteile	194
		7.5.2	Tauschen des Pumpenkopfs	195
		7.5.3	Tauschen der Kolben	200
		7.5.4	Reinigen des Kolbens	202
		7.5.5	Tauschen des Stützrings	203
		7.5.6	Tauschen der Kolbendichtung	208
		7.5.7	Tauschen der Hinterspüldichtung	215
		7.5.8	Tauschen der Pumpenkopfdichtungen	219
		7.5.9	Warten der Pumpenkopf-Kugelventile	224
	7.6	Wartu	ing des Flowmeters	229
		7.6.1	Tauschen des Flowmeter-Filters	229
		7.6.2	Tauschen der Flowmeter-Abfallleitung	231
	7.7	Wartu	Ing der Lösungsmittelschläuche	234
		7.7.1	Entleeren der Lösungsmittelschläuche	234
		7.7.2	Tauschen der Lösungsmittelschläuche vom Behälter zur Pumpe	234
		7.7.3	Tauschen der Lösungsmittelschläuche vom Absperrventil zum Pumpenkopf	239
		7.7.4	Tauschen des Lösungsmittelfilters	240
	7.8	Wartu	Ing der Probenschleife	242
		7.8.1	Komponenten der Probenschleife	242
		7.8.2	Entfernen der Probenschleife	243
		7.8.3	Einbauen der Probenschleife	244
		7.8.4	Abschließen des Probenschleifen-Tauschs	245

7.9	Wartu	ng der Nadeleinheit	247
	7.9.1	Komponenten der Nadeleinheit und Nadelsitz	247
	7.9.2	Sichern der Nadeleinheit	248
	7.9.3	Entsichern der Nadeleinheit	249
	7.9.4	Tauschen des Nadelsitzes und der Nadeleinheit	251
7.10	Wartu	ng des Kopfes der Dosiereinheit	260
	7.10.1	Komponenten des Kopfes der Dosiereinheit	260
	7.10.2	Entfernen und Anschließen von Kapillaren am Kopf der Dosiereinheit.	260
	7.10.3	Tauschen des Kopfes der Dosiereinheit	261
7.11	Wartu	ng der Drainagepumpe	265
	7.11.1	Testen des Drainagepumpen-Schlauchs	265
	7.11.2	Tauschen des Drainagepumpen-Schlauchs	267
7.12	Wartu	ng des Nadel-Waschsystems	270
	7.12.1	Tauschen der Waschschläuche vom Behälter zum Eingang der Nadel-	
		Waschpumpe	270
	7.12.2	Tauschen des Waschschlauchs vom Auslass der Nadel-Waschpumpe zu Waschport	um 272
	7.12.3	Tauschen der Waschschläuche vom Behälter zu den Einlass-Ports am Degaser	273
	7.12.4	Tauschen der Waschschläuche vom Degaser-Ausgang zum Umschaltve	entil 276
	7.12.5	Tauschen des Kugelventils im Nadel-Waschsystem	277
7.13	Wartu	ng der Hinterspülung	279
	7.13.1	Testen der Hinterspülung auf Undichtigkeit	279
	7.13.2	Tauschen des Hinterspülschlauchs zum Behälter	281
	7.13.3	Tauschen des Hinterspül-Einlassschlauchs	284
	7.13.4	Tauschen des Hinterspülpumpen-Einlassschlauchs	287
	7.13.5	Tauschen des Peristaltikschlauchs in der Hinterspülpumpe	289
	7.13.6	Tauschen der Hinterspülschläuche zwischen Hinterspülpumpe und	
		Hinterspüldetektor	292
	7.13.7	Tauschen des Hinterspüldetektors	293
7.14	Prüfen	der Autosampler-Ventile auf Dichtigkeit	296
7.15	Wartu	ng des Säulenthermostaten	298
	7.15.1	Tauschen der Säulen	298
	7.15.2	Visuelle Prüfung der Säulenschaltventile auf Dichtigkeit	299
	7.15.3	Tauschen des Vanquish-Schaltventils	300
	7.15.4	Tauschen des VICI-Schaltventils	303
	7.15.5	Tauschen des Statoren und/oder der Rotordichtung des VICI-Schaltver	ntils 305
	7.15.6	Tauschen des passiven Vorheizers	307

	7.16	Wartu	ng des System Controllers und des Displays	309		
		7.16.1	Tauschen des System Controllers	309		
		7.16.2	Tauschen der SD-Karte	315		
		7.16.3	Tauschen des Displays	317		
	7.17	Tausch	nen des Solvent Racks	319		
	7.18	Tausch	nen des Einsatzes für Flaschen	322		
		7.18.1	Entfernen des Einsatzes für Flaschen	322		
		7.18.2	Einsetzen des Einsatzes für Flaschen	323		
	7.19	Tausch	nen der Sicherungen und Türen	324		
		7.19.1	Tauschen der Sicherungen	324		
		7.19.2	Tauschen der Türen	325		
	7.20	Tausch	nen des Moduleinschubs	327		
		7.20.1	Vorbereiten für das Entfernen der Pumpe	327		
		7.20.2	Entfernen des Moduleinschubs der Pumpe	328		
		7.20.3	Vorbereiten für das Entfernen des Autosamplers	329		
		7.20.4	Entfernen des Moduleinschubs des Autosamplers	331		
		7.20.5	Vorbereiten für das Entfernen des Säulenthermostaten	332		
		7.20.6	Entfernen des Moduleinschubs des Säulenthermostaten	333		
		7.20.7	Zurückschicken des Moduleinschubs	335		
		7.20.8	Installieren des Moduleinschubs	336		
	7.21	Vorbei	reiten des Systems für den Transport	338		
	7.22	Anbrin	gen der Tragegriffe	340		
8	Feh	Fehlersuche				
	8.1	Allgem	neine Informationen zur Fehlersuche	344		
	8.2	Meldu	ngen	346		
		8.2.1	Übergreifende Meldungen	346		
		8.2.2	Meldungen der Pumpe	352		
		8.2.3	Meldungen des Autosamplers	363		
		8.2.4	Meldungen des Säulenthermostaten	368		
	8.3	System	n-Fehlersuche	369		
		8.3.1	Peaks	371		
		8.3.2	Druck	376		
		8.3.3	Retentionszeitschwankungen	377		
		8.3.4	Basislinie	379		
		8.3.5	Längere Ladezeit der Trap-Säule	380		
		8.3.6	Temperaturregelung	380		

		8.3.7	Stromversorgung und Kommunikation (Systemmodule) 381
		8.3.8	Probleme bei System Controller, Display und der Benutzeroberfläche. 383
	8.4	Testen	auf Undichtigkeit
9	Spe	zifikat	tionen
	9.1	Techni	sche Spezifikationen
		9.1.1	Umgebungsbedingungen 388
		9.1.2	Leistungsaufnahme
		9.1.3	Abmessungen
		9.1.4	Emissionsschalldruckpegel 389
		9.1.5	Gewicht
	9.2	Leistur	ngsspezifikationen
		9.2.1	Leistungsspezifikationen des Systems 390
		9.2.2	Leistungsspezifikationen der Pumpe 391
		9.2.3	Leistungsspezifikationen des Autosamplers 392
		9.2.4	Leistungsspezifikationen des System Controllers 394
		9.2.5	Leistungsspezifikationen des Säulenthermostaten 394
10	Zub	ehör,	Ersatzteile und Verbrauchsmaterialien 397
	10.1	Allgem	neine Informationen
	10.2	Zubeh	örkit 399
		10.2.1	System
		10.2.2	Säulenthermostat 402
	10.3	Optior	ales Zubehör 403
	10.4	Ersatzt	teile und Verbrauchsmaterialien 406
		10.4.1	Kapillaren, Kapillarzubehör und Drainage 406
		10.4.2	System-Ersatzteile und Werkzeuge 407
		10.4.3	System Controller-Ersatzteile 408
		10.4.4	Lösungsmittel und Waschsysteme 409
		10.4.5	Pumpen-Ersatzteile 411
		10.4.6	Autosampler-Ersatzteile 413
		10.4.7	Ersatzteile des Säulenthermostaten 414
		10.4.8	Schnittstellen-Kabel und Netzkabel 414

11	Anhang 417	7
	11.1 Informationen zur Konformität	8
	11.1.1 Konformitätserklärungen 41	8
	11.1.2 WEEE-Konformität 41	9
	11.1.3 Einhaltung der FCC-Richtlinien 41	9
	11.1.4 Versionsgeschichte der Anleitung 42	0
	11.2 Digital I/O (Anschlussbelegung) 42	1
	Index	3

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 enthält Anweisungen für Installation, Aufbau, Inbetriebnahme, Außerbetriebnahme, Betrieb, Wartung und Fehlersuche für Ihr Vanquish Neo-System.

Diese Anleitung enthält auch Sicherheitsinformationen, Vorsichtsmaßnahmen und spezielle Hinweise. Halten Sie diese korrekt ein, um Personenschäden, die Beschädigung des Systems oder den Verlust von Daten zu vermeiden.

Hinweise zu den Abbildungen

Die Beschreibungen in dieser Anleitung beziehen sich auf eine Standard-Vanquish™-Neo-Systemkonfiguration.

- Die Abbildungen in dieser Betriebsanleitung dienen dem grundlegenden Verständnis. Sie können vom eigentlichen Modell des Systems oder der Komponente abweichen.
- Aus den Abbildungen in dieser Anleitung können keine Ansprüche hergeleitet werden.

1.2 Konventionen für 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 für einen bestimmten Abschnitt oder eine bestimmte Anweisung gelten, befinden sich in dem jeweiligen Abschnitt oder in der Anweisung, für 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.

1.3 Besondere Hinweise

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 System 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 Systems zu optimieren.

1.4 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

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 auf die Perspektive einer Person, die direkt vor dem System 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)

1.5 Referenzdokumentation

Zusätzlich zu dieser Anleitung stehen weitere Dokumente zur Verfügung.

Thermo Fisher Scientific stellt die aktuellen Betriebsanleitungen als PDF (Portable Document Format) zur Verfügung, Sie können diese von unserer Webseite für Kunden-Dokumentation herunterladen. Um die PDF-Dateien öffnen und lesen zu können, werden Adobe™ Reader™ oder Adobe™ Acrobat™ benötigt.

Gehen Sie zu folgender Webseite: www.thermofisher.com/HPLCmanuals

Hardware-Dokumentation

Informationen zu bewährten Methoden und zur Einrichtung unterstützter Anwendungen Ihres Vanquish Neo-Systems finden Sie im Vanquish Neo System User Guide.

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 Systems.

2.1 Sicherheitssymbole und Signalwörter

2.1.1 Sicherheitssymbole und Signalwörter in dieser Anleitung

Diese Anleitung enthält Sicherheitshinweise zum Schutz von Personen, die das System 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 System installieren oder betreiben, so dass Sie mit dem System und der Anleitung vertraut sind. Die Anleitung enthält wichtige Informationen zur Sicherheit der Anwender sowie zu Gebrauch und Wartung des Systems.
- Bewahren Sie diese Anleitung stets beim System 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 System

In der Tabelle sind die Sicherheitssymbole aufgeführt, die am System oder auf Aufklebern am System angebracht sind. Befolgen Sie die Sicherheitshinweise in dieser Anleitung, um eine Verletzungsgefahr für den Bediener und/oder Schäden am System 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
0	Bezeichnet einen Ein-/Ausschalter, der wechselweise als Ein- oder Aus-Taster für das System wirkt.
(U) VSC	Bezeichnet einen Ein-/Ausschalter, der den System-Controller ein- oder ausschaltet.
\sim	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.
	Weist auf eine mögliche Klemmgefahr hin. Halten Sie Abstand, um Verletzungen an den Händen zu vermeiden.
<u> </u>	Weist auf einen Anschluss mit einer Schutzerdung hin Kann verwendet werden, um eine Vorrichtung zur Schutzerdung anzuschließen, zum Beispiel eine Säule.

2.1.4 Typenschilder

Die Typenschilder sind auf den einzelnen Systemmodulen bei den elektrischen Anschlüssen angebracht. Das Typenschild gibt Auskunft über Seriennummer, Teilenummer, Modulnamen, Revisionsnummer, Netzspannung und Nennleistung der Sicherungen.

Seriennummernaufkleber des Systems

Zusätzlich ist ein Seriennummernaufkleber des Systems unten links in der Schublade des Systemsockels angebracht. Der Seriennummernaufkleber des Systems gibt Auskunft über Seriennummer, Teilenummer, Systemnamen und Revisionsnummer.

Typenschild	
	Ein weiteres Typenschild an der Leckage-Wanne der einzelnen Systemmodule gibt Auskunft über den Modulnamen, die Seriennummer, die Teilenummer und die Revisionsnummer. Für die Kommunikation mit Thermo Fisher Scientific benötigen Sie zur leichteren Identifizierung Ihres Gerätes die Informationen dieses Typenschilds.
System Controller	
	Das Typenschild ist auf der kurzen Seite des System Controllers angebracht. Es gibt Auskunft über Seriennummer und Netzspannung.
Display	
	Das Typenschild ist auf der Rückseite des Displays angebracht. Es gibt Auskunft über Modulnamen, Seriennummer, Teilenummer und Revisionsnummer (wenn vorhanden).

2.2 Verwendungszweck

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

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

Das Vanquish Neo-System und seine Module wurden 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 Neo-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 für 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 Systems die allgemeinen Sicherheitsinformationen in diesem Abschnitt sowie alle anderen in dieser Anleitung aufgeführten spezifischen Sicherheitshinweise und Vorsichtsmaßnahmen beachten.



Wird das System nicht entsprechend den Angaben von Thermo Fisher Scientific eingesetzt, kann dies die im System enthaltenen Schutzvorkehrungen beeinträchtigen. Beachten Sie Folgendes:

- Betreiben Sie das System nur innerhalb der technischen Spezifikationen.
- Verwenden Sie ausschließlich Ersatzteile und zusätzliche Komponenten, Optionen und Peripheriegeräte, die von Thermo Fisher Scientific ausdrücklich für das System autorisiert und freigegeben sind.
- Führen Sie nur die Arbeiten durch, die in dieser Betriebsanleitung und in weiteren Dokumenten für das System beschrieben sind.
 Folgen Sie allen Anweisungen Schritt für Schritt und verwenden Sie die in der Anleitung empfohlenen Werkzeuge.
- Öffnen Sie das Gehäuse des Systems 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 Systems 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 System installieren und/oder bedienen.

Installation



Die Installation des Systems 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 Systems durch eine andere Person als einen Thermo Fisher Scientific-Servicetechniker erfolgen, trägt diejenige Person die Verantwortung dafür, dass die Sicherheit des Systems gewährleistet ist.

Betrieb



Allgemeiner Betrieb

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

Alle Anwender müssen die Gefahren kennen, die von dem System 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

2.3.4 Sicherheitsmaßnahmen beim Umgang mit elektrischen Geräten

Schutzhandschuhe.



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.

2.3.5 Allgemeine Restrisiken

Beachten Sie die folgenden allgemeinen Restrisiken, wenn Sie mit dem System 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.
- Vermeiden Sie den Umgang mit Lösungsmittelbehältern über Kopfhöhe.
- Betreiben Sie das System 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 System nicht in einer Umgebung mit brennbaren Gasen und Dämpfen.



WARNUNG—Brennbare und Gefährliche Dämpfe

Brennbare oder gefährliche Dämpfe können aus schlecht verschlossenen Probengefäßen mit brennbaren oder flüchtigen Proben austreten und sich im Autosampler ansammeln. Dies birgt Gefahren für Gesundheit und Sicherheit und führt zu falschen Ergebnissen.

- Verwenden Sie ausschließlich Probenfläschchen oder Mikrotiterplatten, die mit Hilfe von Verschlusskappen, Verschlussmatten, oder Verschlussfolien gasdicht verschlossen sind. Diese finden Sie in der aktuellen Übersicht der von Thermo Fisher Scientific empfohlenen Verschlüsse.
- Prüfen Sie alle Probenfläschchen vor der Verwendung auf Risse oder Schäden. Verwenden Sie keine Probenfläschchen mit Rissen oder Beschädigungen.



VORSICHT—Austreten von Gefahrstoffen aus PEEK-Kapillaren

Einige Kapillaren im System sind aus PEEK gefertigt. Wenn PEEK-Kapillaren aufquellen oder von Säuren angegriffen werden, können sie undicht werden oder bersten. Bei Kontakt mit einigen Chemikalien, wie zum Beispiel Trichlormethan (CHCl₃), 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 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 die Systemmodule vom Stromnetz.

2.4 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. Die Systemmodule erfüllen bei Auslieferung des Systems alle geltenden Vorschriften hinsichtlich elektromagnetischer Verträglichkeit (EMV) und Sicherheitsstandards.

Änderungen an den Systemmodulen können dazu führen, dass einer oder mehrere dieser EMV- und Sicherheitsstandards nicht mehr eingehalten werden. Änderungen an den Systemmodulen beinhalten auch den Austausch von Teilen oder das Hinzufügen von Komponenten, Zusatzmodulen 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 System hat das Werk in einwandfreiem Zustand verlassen.

3 Übersicht über das System

Dieses Kapitel stellt Ihnen die besonderen Merkmale des Systems und dessen wichtigste Komponenten vor.

3.1 Systemmerkmale

Das System ist wie folgt ausgestattet:

Hauptmerkmale

- Vorinstalliertes System, das vielseitige Möglichkeiten für Workflows bietet
- Betriebsdruck von bis zu 150 MPa
- Flussbereich von 100 nL/min bis 100 μL/min
- Integrierte Funktionen für Autostart und automatisiertes Purgen
- Fortlaufende Systemüberwachung im Hintergrund, um ungeplante Ausfallzeiten zu vermeiden
- Diagnosemöglichkeiten zur schnellen Fehlerbehebung und für verbesserte Verfügbarkeit
- Benutzeroberfläche mit Berührungsbildschirm, über die eine Geräteüberwachung, Steuerung und Wartungsunterstützung möglich ist

Besonderheiten der Pumpe

- Binäre Pumpe mit zwei in Serie geschalteten Kolben je Kanal
- Schwimmend gelagerte Kolben zur Kompensation kleiner mechanischer Toleranzen innerhalb der Spezifikation und damit höhere Robustheit der Pumpe
- Ein Flowmeter f
 ür Flussraten von 100 nL/min bis 100 μL/min; bis auf die Systemkapillaren keine Änderungen an der Hardware erforderlich
- Aktive Hinterspülung für eine hohe Lebensdauer von Kolben und Kolbendichtungen
- Automatisierte Purge-Ventile für einfachen Startup und Diagnose
- Integrierte Funktionen für Autostart und automatisiertes Purgen
- Integrierte automatische Absperrventile verhindern das Austreten von Lösungsmitteln und gewährleisten ein störungsfreies Ansaugen von Lösungsmitteln
- Integrierter Vakuumdegaser für die Nadel-Waschflüssigkeiten des Autosamplers

Besonderheiten des Autosamplers

- Split-Loop-Injektionsprinzip mit geringer Verschleppung und zwei Nadel-Waschflüssigkeiten (schwach und stark)
- Waschen von Nadel, Nadelsitz und Probenschleife von Innen mit schwacher und starker Waschflüssigkeit
- Waschen der äußeren Nadelfläche mit schwacher und starker Waschflüssigkeit für minimale LC/MS-Verschleppung
- Injektionsvolumina von bis zu 100 μL mit nur einem Hub
- Injektionsvolumina über 100 μL durch mehrfaches Aufziehen zum Anreichern größerer Volumina
- Integrierte Funktionen für Autostart und automatisiertes Purgen
- Bodenerkennung, damit kleine Volumina mit minimalem Probenverlust aufgezogen werden können
- Temperaturregelung zur Kühlung bzw. Erwärmung des Probenraums
- Barcodeleser im Autosampler f
 ür die Erkennung von Probenhaltern, Mikrotiterplatten mit Barcodes und des Vanquish-Probenhalterformats
- Aufnahme von Probenhaltern und Mikrotiterplatten mit dem von der Society for Biomolecular Screening definierten Flächenmaß (SBS-Flächenformat)
- Hinterspülung durch den Kopf der Dosiereinheit, um die Lebensdauer der Dichtungen zu erhöhen
- Integrierte Beladungsfunktion

Besonderheiten des Säulenthermostaten

- Säulenraum (34 x 8 x 4 cm, B x H x T) für Säulen, Trap-Säulen und Kapillaren in verschiedenen Workflows
- Platz für bis zu zwei Schaltventilen innerhalb des Säulenraums für fortschrittliche Anwendungen
- Der Säulenraum und die darin befindlichen Komponenten können über einen großen Temperaturbereich thermostatisiert werden; dabei beträgt die minimale Temperatur 5 °C über Umgebungstemperatur und die maximale Temperatur 80 °C ohne Ventile oder 60 °C mit Vanquish Neo-Ventilen mit niedriger Dispersion (LD)

- Optimiertes Design für gleichmäßigen Luftstrom und gleichmäßige Wärmeverteilung
- Schnelle Aufwärmzeit (von 35 °C auf 65 °C ±1 °C in weniger als 12 Minuten)
- Erweiterte Temperaturregelung f
 ür eine Temperaturstabilit
 ät von ±0,05 °C
3.2 Beschreibung des Systems

Das Vanquish-System wurde für die Verwendung in Low-Flow Liquid Chromatography (LC)-Anwendungen und für den Anschluss an ein Massenspektrometer entwickelt.

Das Vanquish-System kann folgende Module enthalten.



Abbildung 1: Systemturm

Nr.	Beschreibung
1	Systemsockel
2	System-Controller
3	Pumpe
4	Autosampler
5	Säulenthermostat (optional)
6	Solvent Rack
7	Display

3.3 Systemsockel

Der Systemsockel trägt die Pumpe, den Autosampler, den System Controller und den optionalen Säulenthermostat.

3.3.1 Komponenten des Systemsockels

Der Systemsockel besitzt:

- Ein-/Ausschalter zum Ein- und Ausschalten aller Systemmodule und des System Controllers
- Schublade zum Aufbewahren von Werkzeug und kleinen Systemkomponenten
- System-Drainage-Ablauf zum Anschluss eines Ablaufschlauchs
- Systemsockelschlüssel zum Umschalten zwischen Verschiebemodus und stationärem Modus



Abbildung 2: Systemsockel (Ansicht von vorne und hinten)

Nr.	Beschreibung
1	Vanquish System Controller Ein- und Ausschalter
	Zum Ein- und Ausschalten des System-Controllers.
2	System-Ein-/Ausschalter Alle Module des Systems, die über einen System-Interlink-Port mit dem Systemsockel verbunden sind, können gemeinsam ein- bzw. ausgeschaltet werden.
3	Schublade für Werkzeug Werkzeug für Installation oder Wartung des Vanquish-Systems kann hier aufbewahrt werden. Die Schublade öffnet und schließt sich mit einem leichten Druck auf die Mitte der Schubladenfront.
4	Reserviert für zukünftige Verwendung
5	System-Drainage-Ablauf Zum Anschluss eines Ablaufschlauchs.

Nr.	Beschreibung
6	Schlösser - auf der linken und rechten Seite des Systemsockels Ermöglicht das Umschalten zwischen Verschiebemodus und stationärem Modus.
7	System-Interlink-Port Ermöglicht den Anschluss der anderen Module zum Ein- und Ausschalten über den Vanquish-Systemsockel.
8	VSC-Port Zum Anschluss eines System-Controllers.

3.3.2 Systemsockelschloss

Der Systemsockel verfügt über einen Sperrmechanismus. Wenn Sie den Systemsockel zum Beispiel bei der Installation oder bei der Wartung verschieben wollen, entsperren Sie ihn zuvor. Um zu verhindern, dass der Systemsockel während des Betriebs auf dem Arbeitstisch verschoben werden kann, sperren Sie den Systemsockel.

Sie können zwischen dem Verschiebemodus und dem stationärem Modus mit den Schlüsseln, die mit dem Sockel geliefert wurden, hinund herschalten.



Abbildung 3: Systemsockelschlüssel

3.4 System Controller

Die Abbildung zeigt die Komponenten des System Controllers.



Abbildung 4: Systemsockel mit VSC-Ein- und Ausschalter und System Controller

Nr.	Beschreibung
1	VSC-Ein- und Ausschalter
2	Federklammern
3	Storage Device (SD) Kartensteckplatz
4	Reserviert für zukünftige Verwendung

3.5 Pumpe

3.5.1 Innenansicht der Pumpe

Die für den Anwender zugänglichen Komponenten der Pumpe befinden sich direkt hinter den Gerätetüren.



Abbildung 5: Innenansicht der Pumpe

Nr.	Beschreibung
1	Tastatur zur Bedienung der Pumpe
2	Pumpenkopf A (linker Pumpenkopf)
3	Clips zur Führung von Lösungsmittelschläuchen zu den Pumpenköpfen
4	Pumpenkopf B (rechter Pumpenkopf)
5	Absperrventile
6	Leaksensor
7	Typenschild mit Angabe von Modulnamen, Seriennummer, Teilenummer und Revisionsnummer
8	Drainage-Ablauf
9	Detektor der Hinterspülung (Tropfendetektor)
10	Degaser für Nadel-Waschflüssigkeiten des Autosamplers
11	Hinterspülpumpe
12	Kanal für die Durchführung der Schläuche, einschließlich Schlauchführungen
13	Flowmeter, mit:
13a	Flowmeter-Einlass B
13b	Pumpenausgang (Flowmeter-Ausgang)
13c	Flowmeter-Abfallleitung (Purge Out)
13d	Flowmeter-Einlass A

2b -_2d __2e -3a A 3f Α -2c ___3b _3c 3d 4 2a 1 3g В -3h B 2 3

3.5.2 Funktionsprinzip der Pumpe



Nr.	Beschreibung
1	Lösungsmittelschläuche: A und B
2	Flussgeregelte Pumpe mit identischen Pumpenblöcken (Lösungsmittelkanälen), A und B
2a	Pumpenblöcke A und B, jeder Pumpenblock mit:
2b	Arbeitszylinder
2c	Arbeitsdrucksensor
2d	Ausgleichszylinder
2e	Primärdrucksensor Der Primärdruck ist der Druck im Pumpenkopf.
3	Flowmeter, je nach Kanal, mit:
3a	Flowmeter-Einlasskapillare (Kapillare vom Pumpenkopf)
3b	Flowmeter-Filter
3c	Automatisches Purge-Ventil mit Waste-Ablauf
3d	Fluss-Sensor
3e	Widerstandskapillare
3f	Systemdrucksensor
3g	Flowmeter-Ausgang mit fluidischem T-Stück (Pumpenausgang)
3h	Waste-Auslass
4	Kapillare zum System

Die Pumpe ist eine binäre Pumpe mit Flowmeter zur Flussregelung:

- Jeder Pumpenblock hat einen Arbeitszylinder und einen Ausgleichszylinder. Die Zylinder sind in Serie geschaltet. Das Lösungsmittel durchläuft die Zylinder nacheinander.
- Jeder Pumpenblock liefert den Fluss kontinuierlich.
 Der Arbeitszylinder fördert, während der Ausgleichszylinder gleichzeitig gefüllt wird. Der Ausgleichszylinder dient als Speicher und fördert, während der Arbeitszylinder den Saughub ausführt.
- Die Sensoren im Flowmeter messen und regeln die Teilflüsse der Pumpenblöcke.
- Die Teilflüsse werden im fluidischen T-Stück am Flowmeter-Ausgang zusammengeführt. Der zusammengeführte Fluss verlässt das Flowmeter mit dem eingestellten Gesamtfluss und der vorgegebenen Lösungsmittelzusammensetzung.

3.6 Autosampler

3.6.1 Innenansicht des Autosamplers

Die für den Anwender zugänglichen Komponenten des Autosamplers befinden sich direkt hinter den Gerätetüren.



Abbildung 7: Innenansicht VN-A10 Autosampler

Nr.	Beschreibung
1	Probenraum mit Drehteller
2	Drainagepumpe
3	Leckage-Wanne mit:
3a	Tastatur mit Statusanzeigen
3b	Führungsloch für die Kapillare zur Pumpe
3c	Typenschild mit Angabe von Modulnamen, Seriennummer, Teilenummer und Revisionsnummer
3d	Kanal zum Drainage-System
3e	Drainage-Abläufe
3f	Leaksensor
4	Nadel-Waschpumpe
5	Injektionseinheit (Die Detailansicht finden Sie auf der nächsten Seite)
6	Isolierabdeckung
7	Einlass-Port der Hinterspülung
8	Führungen für die Kapillare zum Säulenthermostat



Injektionseinheit des Autosamplers

Abbildung 8: Detailansicht der Injektionseinheit im VN-A10 ohne Isolierabdeckung

Nr.	Beschreibung
1	Transportsicherungsschraube für die Nadeleinheit
2	Nadeleinheit mit Niederhalter
3	Probenschleifeneinheit
4	Nadelsitz
5	Injektionsventil
6	Lösungsmittelventil
7	Kugelventil
8	Umschaltventil
9	Kopf der Dosiereinheit
10	Drucksensor

3.6.2 Funktionsprinzip des Autosamplers

Das Funktionsprinzip des Autosamplers basiert auf dem Split-Loop -Injektionsprinzip. Bei diesem Injektionsprinzip ist die Nadel Teil der Probenschleife. Im Folgenden wird der Aufbau für direkte Injektion beschrieben.

Aufziehen der Probe, Ventile in Bypass-Position

- Zu Beginn des Injektionsvorgangs f\u00e4hrt die Probennadel vom Nadelsitz zur vordefinierten Position der Probe auf dem Drehteller. Dort taucht die Nadel in das Probengef\u00e4\u00df ein.
- Die Dosiereinheit zieht die Probe vom Probengefäß durch die Probennadel in die Probenschleife. Anschließend fährt die Probennadel vom Probengefäß hoch.
- Um zwischen den Injektionen Probenreste von der Probennadel zu spülen, kann das Äußere der Nadel gewaschen werden. In diesem Fall fährt die Probennadel in den Waschport und taucht ein. Die Nadel-Waschpumpen fördern Waschflüssigkeit zum Waschport, um die Probennadel von außen zu waschen; die Waschflüssigkeit geht dabei in den Waschabfall.
- Die Probennadel fährt zurück zum Nadelsitz, der die Nadelspitze gegen den Systemdruck abdichtet.

Injektion der Probe, Ventile in Inject-Position (siehe Abbildung)

 Das Injektionsventil schaltet auf die Inject-Position, um den benötigten Flussweg herzustellen, sodass das Lösungsmittel aus der Pumpe durch die Probenschleife, den Nadelsitz und die Probennadel fließt und die Probe durch das Injektionsventil auf die Säule spült.

Beginn des Gradienten, Ventile in Bypass-Position

 Wenn die Probe vollständig aus dem Autosampler herausgedrückt wurde, schaltet das Injektionsventil zurück in die Bypass-Position. Der Gradient kann nun mit dem kleinstmöglichen Gradientenverzögerungsvolumen gefördert werden.
 ANMERKUNG: Falls gewünscht kann der Autosampler so konfiguriert werden, dass das Injektionsventil in der Inject-Position bleibt und der Gradient sofort gestartet wird. Da die Probenschleife vollständig mit einbezogen ist, erhöht sich das Gradientenverzögerungsvolumen.

Waschvorgang, Ventile in Wasch-Position

Während das Injektionsventil in Bypass-Position steht, schaltet das Lösungsmittelventil in die Wasch-Position und die Probennadel fährt vom Nadelsitz in den Waschport und taucht ein. In dieser Position werden das Nadeläußere und das Nadelinnere mit starker Nadel-Waschflüssigkeit aus dem Waschport (7) und schwachem Nadel-Waschlösungsmittel aus dem Online-Degaser in der Pumpe (13) gewaschen.



Abbildung 9: Funktionsprinzip VN-Autosampler, Inject-Position

Nr.	Beschreibung
1	Drehteller
2	Probenschleife
3	Nadel
4	Nadelsitz
5	Waschport
6	Nadel-Waschpumpe
7	Nadel-Waschflüssigkeiten (schwach und stark)
8	Injektionsventil
9	Lösungsmittelventil
10	Waschabfall
11	Drucksensor
12	Dosiereinheit
13	Wasch-Lösungsmittel vom Degaser in der Pumpe (schwach und stark)
14	Umschaltventil
15	Pumpe
16	Säule

3.6.3 Anschlussbelegung der Autosampler-Ventile

Im Aufbau für direkte Injektion sind die Ports des Injektionsventils folgendermaßen belegt:



Abbildung 10: Injektionsventil mit angeschlossenen Kapillaren

Port	Angeschlossene Komponente
1	Viper-Verschlussfitting
2	Lösungsmittelventil
3	Nadelsitz
4	Trennsäule (falls ein Säulenthermostat installiert ist)
5	Probenschleife
6	Dosiereinheit
7	Pumpe



Für die direkte Injektion und Workflows vom Typ Trap und Elution sind die Ports des Lösungsmittelventils folgendermaßen belegt:



Port	Angeschlossene Komponente
1	Viper-Verschlussfitting
2	Abfall
3	Injektionsventil
4	Viper-Verschlussfitting
5	Kugelventil
6	Drucksensor
7	Abfall

3.6.4 Probenraum

Der Probenraum des Autosamplers umfasst einen Drehteller mit vier farblich gekennzeichneten Segmenten; wobei jedes Segment zusätzlich mit einem Probenfläschchen-Halter für drei 10 mL-Probenfläschchen ausgestattet ist.

Der Autosampler kann Probenhalter und Mikrotiterplatten mit dem von der Society of Biomolecular Screening definierten Flächenmaß (SBS-Flächenmaß) aufnehmen.

Der Probenraum verfügt über folgende Ausstattungsmerkmale:

- Probenraum-Thermostatisierung, siehe Probenraum-Thermostatisierung (> Seite 50).
- Erkennung und Verifizierung des Probenhalterformats, Erkennung von leeren Segmenten und Bestandsmanagement, siehe Probenhalterformat-Erkennung (> Seite 51).

3.6.4.1 Probenraum-Thermostatisierung

Die Temperaturregelung kann zur Probenraum-Thermostatisierung eingeschaltet werden. Zirkulierende Luft kühlt bzw. erwärmt den Probenraum auf die eingestellte Temperatur und ermöglicht dadurch eine präzise Temperierung der Proben.

Um eine optimale Thermostatisierung zu erzielen, halten Sie die Autosamplertüren während der Probenanalyse geschlossen. Öffnen Sie die Gerätetüren nur, wenn erforderlich, zum Beispiel, um Probengefäße in den Autosampler zu stellen.

Die Probenraum-Thermostatisierung verfügt über einen Temperaturbereich von +4 °C bis +40 °C. Das Kühlen der Proben ist bis maximal 23 K unterhalb der Umgebungstemperatur möglich.

Abhängig von der Umgebungsfeuchtigkeit kann Kondensation während der Probenkühlung auftreten. Eine Drainagepumpe (Peristaltikpumpe) im Autosampler entfernt jede entstehende Kondensationsflüssigkeit aus dem Probenraum.

Der Autosampler ist mit einer Isolierabdeckung ausgestattet, die die Probenschleife abdeckt. Diese Abdeckung schützt den Innenraum vor hoher Feuchtigkeit und übermäßiger Kondensation, wenn der Probenraum thermostatisiert wird. Wenn die Abdeckung nicht ordnungsgemäß angebracht ist, kann die Thermostatisierung nicht eingeschaltet werden bzw. wird automatisch abgeschaltet.

3.6.4.2 Drehteller

Der Drehteller ist in vier farblich gekennzeichnete Segmente eingeteilt: rot (**R**), grün (**G**), blau (**B**) und gelb (**Y**).

Jedes Segment ist in folgende Bereiche aufgeteilt:

- Ablagefläche für einen Probenhalter oder eine Mikrotiterplatte
- Zusatz-Fläschchenhalter mit drei Positionen für Probenfläschchen mit einem Volumen von bis zu 10 mL, wie Reagenzfläschchen Der Zusatz-Fläschchenhalter befindet sich in den jeweiligen Segmenten links von der Ablagefläche für Probenhalter bzw. Mikrotiterplatten.



Abbildung 12: Detailansicht eines Segments im Drehteller

Nr.	Beschreibung
1	Barcode-Label 'leer'
	Label, das dem Barcodeleser anzeigt, dass das Segment leer ist
2	Farbkennzeichnung für das Segment (hier: R für das rote Segment)
3	Ablagefläche im Segment Fläche, um einen Probenhalter oder eine Mikrotiterplatte zu platzieren Mit Ausrichtungs-Ecken und -Punkten, um den Probenhalter oder die Mikrotiterplatte auszurichten
4	Zusatz-Fläschchenhalter Für drei Probenfläschchen mit einem Volumen von bis zu 10 mL
5	Farbkennzeichnung für den Zusatz-Fläschchenhalter (<i>hier:</i> R1 zum Beispiel für die erste Fläschchenhalter-Position im roten Segment)

3.6.4.3 Probenhalterformat-Erkennung

Ein Barcodeleser im Probenraum ermöglicht die automatische Erkennung von Probenhaltern und Mikrotiterplatten, die mit einem speziellen Vanquish-2D-Barcode zur Format-Erkennung bedruckt sind.

Während des Betriebs führt der Barcodeleser eine Bestandsaufnahme (engl. 'Inventory Scan') durch und liest die Vanquish-Probenhalterformat-Barcodes. Der Barcodeleser identifiziert dabei automatisch das Probenhalter-Format und die Ausrichtung. Diese Information wird an das Chromatographie-Datensystem gesendet. Um die Erkennung des Probenhalter- und Mikrotiterplatten-Formats mit dem Barcodeleser im Autosampler zu ermöglichen, verwenden Sie Probenhalter und Mikrotiterplatten, die mit einem solchen Barcode zur Format-Erkennung bedruckt sind.

In jedem Segment befindet sich ein Barcode-Aufkleber, der dem Barcodeleser anzeigt, dass das Segment leer ist, wenn weder ein Probenhalter noch eine Mikrotiterplatte eingesetzt ist.



Abbildung 13: Probenhalter mit Barcodes zur Format-Erkennung

Nr.	Beschreibung
1	Probenposition A1
2	Vanquish-Probenhalterformat-Barcode

3.7 Säulenthermostat

3.7.1 Innenansicht des Säulenthermostaten

Die für den Anwender zugänglichen Komponenten des Säulenthermostaten befinden sich direkt hinter den Gerätetüren.



Abbildung 14: Innenansicht des Säulenthermostaten

Nr.	Beschreibung
1	Tastatur
2	Linkes Säulenschaltventil (falls installiert; hier: Vanquish-Ventil)
3	Säulen-ID-Ports
4	Lüfter
5	Temperatursensor
6	Kapillarführungen
7	Rechtes Säulenschaltventil (falls installiert; hier: Vanquish-Ventil)
8	Typenschild mit Angabe von Modulnamen, Seriennummer, Teilenummer und Revisionsnummer
9	Einbaupositionen für den passiven Vorheizer
10	Säulenhaltespangen

3.7.2 Funktionsprinzip des Säulenthermostaten

Wichtigste Anforderung an einen Säulenthermostaten für die Flüssigkeitschromatographie ist, dass die eingestellte Temperatur so konstant wie möglich gehalten wird. Durch das moderne Design des Säulenthermostaten ist es möglich, die Temperatur im Säulenraum auf ±0,1 °C konstant zu halten.

Der Temperaturstabilität kommt für das Erzielen guter Messergebnisse dabei eine größere Bedeutung zu, als der Temperaturgenauigkeit. Der Säulenthermostat bietet eine Temperaturstabilität von ±0,05 °C.

Heizelemente erwärmen die Luft im Säulenraum und die sich darin befindenden Komponenten auf die eingestellte Temperatur. Das optimierte Design ermöglicht einen gleichmäßigen Luftstrom.

Dadurch wird Folgendes sichergestellt:

- Die Temperatur der stationären Phase bleibt über die gesamte Säulenlänge gleich.
- Die Säule und der Eluent besitzen während der Trennung dieselbe Temperatur.
- Die analytische Trennung erfolgt bei der eingestellten Temperatur.

3.7.3 Übersicht der Säulenschaltventile

Der Säulenthermostat kann mit einem oder zwei Schaltventilen für erweiterte Anwendungen, einschließlich Säulenschaltung, ausgestattet sein. Sie können die Ventile entsprechend Innenansicht des Säulenthermostaten (> Seite 53) installieren.

Die folgenden Schaltventiltypen können installiert werden:

- Vanquish-Schaltventile
- VICI[™]-Schaltventile

Eine gemischte Konfiguration aus Vanquish-Schaltventilen und VICI-Schaltventilen ist möglich. Sie können Vanquish-Schaltventile an jeder Position installieren, an der kein VICI-Ventilantrieb installiert ist.

Wenn kein Ventil installiert ist, verschließt eine Abdeckung die Öffnung.

Vanquish-Schaltventile

Vanquish-Schaltventile bestehen aus einem Teil. Der Ventilantrieb ist in die Ventileinheit integriert.



Abbildung 15: Vanquish-Schaltventil: Vorderansicht (links), Rückansicht (rechts)

Für die Installation eines Vanquish-Ventils, siehe Installieren des Vanquish-Schaltventils (> Seite 88).

Für den Tausch eines Vanquish-Ventils, siehe Tauschen des Vanquish-Schaltventils (▶ Seite 300).

VICI-Schaltventile

Das VICI-Ventil besteht aus zwei Teilen, dem Ventilkopf (Nr. 1) und dem Ventilklemmring (Nr. 2). Der Ventilklemmring verbindet den Ventilkopf mit dem Ventilflansch (Nr. 3) des Ventilantriebs. Der Ventilantrieb ist nicht in die Ventileinheit integriert. Der Ventilantrieb befindet sich im Gehäuse und ist nicht für den Kunden zugänglich.

Sie können nur das VICI-Ventil installieren oder tauschen. Um den VICI-Ventilantrieb zu installieren, wenden Sie sich an den Kundendienst von Thermo Fisher Scientific.



Abbildung 16: Übersicht des VICI-Ventils (links); Ventilöffnung mit Flansch (rechts)

Nr.	Beschreibung
1	Kopf
2	Klemmring
3	Flansch (im Gehäuse)

Für die Installation eines VICI-Ventils, siehe Installieren des VICI-Schaltventils (Seite 90).

Für den Tausch eines VICI-Ventils, siehe Tauschen des VICI-Schaltventils (
 Seite 303).

3.8 Solvent Rack



Die Abbildung zeigt die Komponenten des Solvent Racks.

Abbildung 17: Solvent Rack-Komponenten

Nr.	Beschreibung	
1	Solvent Rack	
1a	Magnetische Vorderseite An der Vorderseite des Solvent Racks können magnetische Etiketthalter angebracht werden, zum Beispiel, um das System mit einem kundenspezifischen Systemnamen zu kennzeichnen. Etiketten sind im Zubehörkit enthalten.	
2	Einsatz für Flaschen ¹	
2a	5 Mulden für 1-L-Behälter	
2b	2 Mulden für 2-L-Behälter	
2c	4 Mulden für 0,25-L-Behälter	
¹ : Der Einsatz für Flaschen lässt sich leicht herausnehmen. Wenn Sie den Einsatz für		

Flaschen herausnehmen, verwenden Sie vorzugsweise nur die Flüssigkeitsbehälter, die mit dem System mitgeliefert werden. Wenn Sie andere Flüssigkeitsbehälter verwenden, zum Beispiel mit einer größeren Kapazität, stellen Sie sicher, dass der Systemturm stabil steht und nicht kippt. Wenn Sie Zweifel an der Kippsicherheit haben, installieren Sie eines der Kits zur Stabilisierung (siehe Zubehörkit (\triangleright Seite 399)).

3.9 Leakerkennung

Da Undichtigkeiten ein potentielles Sicherheitsrisiko darstellen, sind die Pumpe und der Autosampler mit einem Leaksensor ausgestattet.

Die Leaksensoren im System überwachen die fluidischen Verbindungen auf Undichtigkeiten. Die Flüssigkeit sammelt sich in den Leckage-Wannen und wird zum Ablauf geleitet. Durch den Ablauf fließt die Flüssigkeit über das Drainagesystem in den Abfall.

Wenn ein Leaksensor auf eine Undichtigkeit anspricht, leuchtet die Statusanzeige des betreffenden Moduls rot und ein akustisches Signal ertönt.

Ist eine Undichtigkeit in der Pumpe nicht nach 3 Minuten behoben, stoppt der Pumpenfluss.

Sehen Sie dazu auch

Testen auf Undichtigkeit (Seite 385)

3.10 Nadel-Waschsystem (Übersicht)

Mit Hilfe des Nadel-Waschsystems im Autosampler kann die Nadel von Innen und Außen gewaschen werden, um Probenreste von der Probennadel zu spülen. Die Nadel-Waschzyklen können Teil der Probensequenz sein oder manuell durchgeführt werden.

Die Komponenten des Nadel-Waschsystems innerhalb des Autosamplers sind bei Auslieferung vorinstalliert.

Waschen des Nadeläußeren

Das Nadel-Waschsystem für das Waschen des Nadeläußeren besitzt zwei Waschpumpen, die eine starke oder eine schwache Nadel-Waschflüssigkeit fördern können.

Das Waschen des Nadeläußeren kann automatisch nach dem Aufziehen der Probe oder manuell durchgeführt werden.

Waschen des Nadelinneren

Beim Waschen des Nadelinneren wird das Nadeläußere ebenfalls mitgewaschen. Um die Verschleppung niedrig zu halten, werden auch der Nadelsitz und die Probenschleife mit gewaschen. Zusätzlich zu den beiden Nadel-Waschpumpen für das Waschen des Nadeläußeren wird für das Waschen des Nadelinneren eine schwache Nadel-Waschflüssigkeit verwendet, die durch den Online-Degaser in der Pumpe geleitet wird.



Abbildung 18: Komponenten des Nadel-Waschsystems

Nr.	Beschreibung	
1	Waschschlauch mit Beschriftung SWP (rot) vom Behälter für die starke Nadel-Waschflüssigkeit zum Einlass der Nadel-Waschpumpe, zum Waschen des Nadeläußeren und -inneren mit starker Waschflüssigkeit.	
	Waschschlauch mit Beschriftung WWP (grün) vom Behälter für die schwache Nadel-Waschflüssigkeit zum Einlass der Nadel-Waschpumpe, zum Waschen des Nadeläußeren und -inneren mit schwacher Waschflüssigkeit.	
2	Waschschlauch mit Beschriftung S (rot) vom Behälter für die starke Nadel- Waschflüssigkeit zum Degaser-Eingang in der Pumpe, zum Purgen und für Prozeduren zur kurzzeitigen Außerbetriebnahme.	
	Waschschlauch mit Beschriftung W (grün) vom Behälter für die schwache Nadel-Waschflüssigkeit zum Degaser-Eingang in der Pumpe, zum Waschen des Nadelinneren und zum Purgen.	
3	Waschport	
4	Waschschläuche vom Auslass der Nadel-Waschpumpe zum Waschport	
5	Nadel-Waschpumpe	
6	Waschschläuche vom Degaser-Ausgang an der Pumpe zum Autosampler.	
7	Ablaufschlauch des Nadel-Waschsystems	

3.11 Hinterspülung (Übersicht)

Die Hinterspülung trägt dazu bei, die Lebensdauer von Kolben und Kolbendichtungen zu verlängern; ein Auskristallisieren und die Ablagerung von Partikeln auf den Kolben und Kolbendichtungen werden verhindert, wodurch andernfalls die Komponenten beschädigt werden könnten. Der Flussweg durch das Hinterspülsystem führt durch den Kopf der Dosiereinheit im Autosampler und die Pumpenköpfe in der Pumpe. Die Hinterspülung besteht aus den folgenden Teilen:

- In der Pumpe: Hinterspülpumpe (Peristaltikpumpe), Hinterspülschläuche und Hinterspüldetektor (Tropfendetektor)
- Im Autosampler: Hinterspülschläuche
- Behälter für die Hinterspülflüssigkeit

Bei Auslieferung sind im System die Komponenten der Hinterspülung und die Schlauchverbindungen installiert.



Abbildung 19: Übersicht Hinterspülung

Nr.	Beschreibung
1	Einlass-Port der Hinterspülung (rechter Port)
2	Hinterspülzufuhrleitung vom Behälter zum Einlass-Port

Nr.	Beschreibung
3	Hinterspül-Einlassschlauch vom Einlass-Port zum Kopf der Dosiereinheit
4	Kopf der Dosiereinheit
5 + 6	Hinterspülpumpen-Einlassschlauch
7	Hinterspülpumpe (Peristaltikpumpe)
8	 Hinterspülschläuche in der Pumpe (dünne Schläuche), mit Hinterspülschlauch von der Hinterspülpumpe zum linken Pumpenkopf Hinterspülschlauch am linken Pumpenkopf Hinterspülschlauch zwischen den Pumpenköpfen Hinterspülschlauch am rechten Pumpenkopf Hinterspül-Auslassschlauch von rechtem Pumpenkopf zum Hinterspüldetektor
9	Detektor der Hinterspülung (Tropfendetektor)

Sehen Sie dazu auch

Füllen des Behälters für die Hinterspülflüssigkeit () Seite 133)

3.12 Ein- und Ausschalten

Vorhandene Ein-/Ausschalter und Hauptnetzschalter

Die Abbildung zeigt die Ein-/Ausschaltelemente des Systems.



Abbildung 20: Ein-/Ausschalter und Hauptnetzschalter

Nr.	Beschreibung
1	System-Ein-/Ausschalter
2	Ein- und Ausschalter des System-Controllers
3	Hauptnetzschalter

System-Ein-/Ausschalter

Für eine einfachere Bedienung können Sie den System-Ein-/Ausschalter zum gleichzeitigen Ein-/Ausschalten aller Systemmodule (mit Ausnahme des System-Controllers) verwenden, wenn die Hauptnetzschalter der Systemmodule eingeschaltet sind. Beachten Sie folgende Hinweise:

Status	Beschreibung
Ein	 Alle Systemmodule des Systems, die über einen System-Interlink-Port mit dem Systemsockel verbunden sind, werden durch Drücken des System-Ein/Ausschalters gleichzeitig eingeschaltet.
	 Sie können ein Systemmodul nicht über den System-Ein- /Ausschalter einschalten, wenn der Hauptnetzschalter des Systemmoduls ausgeschaltet ist.
Aus	 Alle Systemmodule des Systems, die über einen System-Interlink-Port mit dem Systemsockel verbunden sind, werden durch Drücken des System-Ein/Ausschalters gleichzeitig ausgeschaltet.
	• Das Drücken des System-Ein-/Ausschalters reicht nicht aus, um die Systemmodule vollständig auszuschalten. Um ein Systemmodul vollständig auszuschalten, müssen Sie es über den Hauptnetzschalter des Systemmoduls ausschalten. Schalten Sie ein Systemmodul über den Hauptnetzschalter aus, wenn Sie dazu aufgefordert werden, zum Beispiel für Wartungsarbeiten.

Ein- und Ausschalter des System-Controllers

Der Ein-/Ausschalter des System-Controllers ist komplett unabhängig vom System-Ein-/Ausschalter. Wenn Sie den System-Controller einschalten, schaltet sich das Display automatisch ein. Nach einem Stromausfall schaltet sich der System-Controller selbständig wieder ein.

Um den System-Controller und das Display (wenn vorhanden) auszuschalten, schalten Sie den Ein-/Ausschalter des System-Controllers aus und warten Sie, bis die Status-LED (Light Emitting Diode) erlischt.

3.13 Betrieb des Systems

Wartung und Systemüberwachung über das Thermo Scientific Vanquish User Interface

Über das Thermo Scientific Vanquish User Interface können Sie das System einrichten und überwachen. Über diese Oberfläche können Sie schnell den Zustand des Systems prüfen, Wartungsprozeduren durchführen, und Probleme lösen. Dies alles ist mit oder ohne Verbindung zum Chromatographie-Datensystem möglich.

Alle Beschreibungen in dieser Bedienungsanleitung zum Thema Software beziehen sich, wenn nicht anders angegeben, auf das Vanquish User Interface.

Einen Überblick über das Vanquish User Interface finden Sie unter Vanquish User Interface (▶ Seite 155).

Betrieb mit einem Chromatographie-Datensystem

Weitere Fragen zum Anschluss und zur Steuerung des Systems an ein Chromatographie-Datensystem beantwortet Ihnen gerne der Thermo Fisher Scientific Kundendienst.

3.14 Purgen (Übersicht)

Purging beschreibt das Spülen des Systems oder von Systemkomponenten für eine bestimmte Zeit mit einer höheren Flussrate. Erforderlich ist Purgen in verschiedenen Situationen und aus unterschiedlichen Gründen. Die Situation und der Grund bestimmen, welches Skript jeweils das beste ist. In den Tabellen finden Sie Informationen zu den häufigsten Situationen. Sie starten die Skripte vom Vanquish User Interface. Folgen Sie gegebenenfalls den Anweisungen des Skripts.

Während der Installation, vor dem Betrieb

Grund	Skript	Zusätzlicher Parameter
Bei der Vorbereitung des Systems für den ersten Betrieb	A02 - Autostart	Keine
Zum Neustart des Systems nach einer langfristigen Außerbetriebnahme	A02 - Autostart	Keine

Während des Betriebs

Grund	Skript	Zusätzlicher Parameter
Zum Austausch eines Lösungsmittelbehälters oder eines Behälters mit der Nadel- Waschflüssigkeit des Autosamplers, wenn der Behälter nicht mehr genug Flüssigkeit für die nächste Analyse enthält (gleicher Typ Flüssigkeit)	B01 - Flüssigkeiten/ Lösungsmittel ändern	Wählen Sie das betroffene Lösungsmittel oder die Nadel-Waschflüssigkeit aus und wählen Sie Nur aktualisieren = An.
Eine oder mehrere Lösungsmittelschläuche oder Schläuche für die Nadel-Waschflüssigkeit des Autosamplers waren leer	B01 - Flüssigkeiten/ Lösungsmittel ändern	Wählen Sie das betroffene Lösungsmittel oder die Nadel-Waschflüssigkeit aus und wählen Sie Nur aktualisieren = Aus.
Zur Umstellung auf einen anderen Typ Flüssigkeiten (Lösungsmittel oder Nadel-Waschflüssigkeit für den Autosampler), zum Beispiel bei Umstellung auf eine andere Analysenmethode	B01 - Flüssigkeiten/ Lösungsmittel ändern	Wählen Sie das betroffene Lösungsmittel oder die Nadel-Waschflüssigkeit aus und wählen Sie Nur aktualisieren = Aus.

Während der Wartung

Grund	Skript	Zusätzlicher Parameter
Pumpenwartung: Zu Beginn einer Wartungsprozedur: Um schädliche Substanzen auszuspülen	CO2 - Pumpe purgen	Was gepurgt werden soll = Gesamte Pumpe Intensität = Standard
Am Ende einer Wartungsprozedur: Um Luftblasen zu entfernen, die sich gegebenenfalls durch die Wartungsprozedur im System befinden		
Autosamplerwartung: Um Luftblasen zu entfernen, die sich gegebenenfalls durch die Wartungsprozedur im System befinden	C04 - Sampler purgen	Was gepurgt werden soll = Gesamter Sampler Intensität = Intensiv

Während der Fehlersuche

Warum	Skript
Um Luftblasen zu entfernen, die sich gegebenenfalls im System befinden, und zum Beispiel zu Pulsation während des Betriebs, starkem Rauschen oder einer nicht-reproduzierbaren Analyse führen	Abhängig von der möglichen Ursache. Siehe System-Fehlersuche (▶ Seite 369).

4 Auspacken

In diesem Kapitel finden Sie Informationen zum Auspacken des großen Versandkartons mit dem vorinstallierten Basissystemturm. Informationen zum Auspacken des optionalen Säulenthermostaten finden Sie unter Auspacken des Säulenthermostaten (▶ Seite 72). Außerdem finden Sie Informationen zum Lieferumfang.

4.1 Beschädigte Verpackung, Mängel bei Ankunft des Geräts

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

Besteht der Verdacht, dass der Inhalt 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.

4.2 Lieferumfang

Im Lieferumfang enthalten sind:

- Vorinstalliertes System
- Display
- Betriebsanleitung des Systems (kann von der Webseite für Kunden-Dokumentation heruntergeladen werden)
- Netzkabel
- Zubehörkit Informationen zum Inhalt des Kits finden Sie in Zubehörkit (>> Seite 399).

4.3 Auspacken des Systems



VORSICHT—Schweres und unhandliches System

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

- Um das System zu greifen und zu transportieren, sind zwei Personen erforderlich, zum Beispiel, um es anzuheben oder zu verschieben.
- Verwenden Sie für den Transport des Systems die Tragegriffe, die am System montiert sind. Transportieren Sie oder heben Sie das System niemals an den Gerätetüren. Dies führt zu Beschädigungen an den Türen oder an den Systemmodulen. Überprüfen Sie, dass die Tragegriffe korrekt angeschraubt sind und nicht herausrutschen können.

ACHTUNG

Selbst wenn Sie den Vanquish-Systemturm nur geringfügig auf dem Arbeitstisch verschieben, kann der Systemsockel beschädigt und dessen Funktionalität beeinträchtigt werden.

Bevor Sie den Systemturm verschieben, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln.

Erforderliche Werkzeuge

Schraubendreher, Torx[™] T30

Gehen Sie wie folgt vor

- 1. Transportieren Sie das System zum Aufstellungsort.
- 2. Entfernen Sie alle Umreifungsbänder vom Karton.
- 3. Heben Sie den Deckel an und entfernen Sie ihn.
- 4. Entnehmen Sie das Zubehör.
- 5. Entnehmen Sie das obere Schaumstoffteil.
- 6. Heben Sie das System aus dem Karton:
 - a) Greifen Sie das System an den Tragegriffen.
 - b) Heben Sie das System langsam und vorsichtig aus dem Versandkarton heraus.
- 7. Stellen Sie das System auf eine stabile Oberfläche und lassen Sie etwaige Schutzfolien noch auf den Oberflächen kleben.

- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (> Seite 73).
- 9. Entfernen Sie sonstiges Verpackungsmaterial.
- 10. Optional: Entfernen Sie die Tragegriffe:
 - a) Lösen Sie die Befestigungsschrauben an den Tragegriffen, bis Sie die Tragegriffe in der Schiene bewegen können. Entfernen Sie die Schrauben nicht vollständig von den Tragegriffen.
 - b) Ziehen Sie die Tragegriffe zur Rückseite des Systems hin aus den Schienen heraus.





11. Ziehen Sie die Schrauben der Tragegriffe an.

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

12. Entfernen Sie die Schutzfolien soweit vorhanden von allen Oberflächen.

4.4 Auspacken des Säulenthermostaten

Beschädigte Verpackung, Mängel bei Ankunft des Geräts

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

Besteht der Verdacht, dass der Inhalt 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

ACHTUNG

Greifen Sie das Gerät zum Anheben oder Verschieben an den Tragegriffen. Transportieren oder heben Sie das Gerät niemals an den Gerätetüren. Dies führt zu Beschädigungen an den Türen oder am Gerät.

Gehen Sie wie folgt vor

- 1. Stellen Sie den Versandkarton auf den Boden und öffnen Sie ihn.
- 2. Entnehmen Sie das Zubehör.
- 3. 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.
- 4. Stellen Sie das Gerät auf eine stabile Oberfläche.
- Entfernen Sie sonstiges Verpackungsmaterial. Einige Oberflächen wie die Türen des Geräts sind für den Transport durch Folien geschützt. 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 Einbau des Säulenthermostaten in den Systemturm (▶ Seite 86).
- 7. Entfernen Sie die Schutzfolien soweit vorhanden von allen Oberflächen.
4.5 Entsperren des Systemsockels

1. Stecken Sie einen Schlüssel in das Schloss auf der einen Seite des Systemsockels.



Abbildung 22: Einstecken der Systemsockelschlüssel

- 2. Stecken Sie einen Schlüssel in das Schloss auf der anderen Seite des Systemsockels.
- 3. Entsperren Sie den Systemsockel, indem Sie die Schlüssel wie unten gezeigt drehen.

In der entsperrten Position können die Schlüssel nicht aus dem Systemsockel gezogen werden.



Abbildung 23: Entsperren des Systemsockels

- 4. Verschieben Sie das System.
- Sperren Sie den Systemsockel (siehe Sperren des Systemsockels (> Seite 118)).

5 Installation

In diesem Kapitel finden Sie Informationen zu den Anforderungen an den Aufstellungsort sowie zum Aufbau, zur Installation und zur Konfiguration des Systems.

5.1 Sicherheitshinweise für die Installation

Beachten Sie folgende Sicherheitshinweise:



Beachten Sie alle Warnhinweise und Vorsichtsmaßnahmen in Sicherheitsmaßnahmen (▶ Seite 26).



WARNUNG—Scharfe Spitze der Injektionsnadel

Die Injektionsnadel besitzt eine scharfe Spitze, die Verletzungen der Haut verursachen kann.

Berühren Sie niemals die Nadelspitze, um Personenschäden zu vermeiden.



VORSICHT—Schweres und unhandliches Gerät

Die System-Module sind zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden an den System-Modulen zu vermeiden, beachten Sie folgende Hinweise:

- Um die Module zu greifen und zu transportieren, sind zwei Personen erforderlich, zum Beispiel, um sie anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn die Module in den Systemturm gehoben oder aus dem Systemturm entfernt werden.
- Verwenden Sie für den Transport der Module die Tragegriffe, die mit den Modulen mitgeliefert wurden. Transportieren Sie oder heben Sie die Module niemals an den Gerätetüren. Dies führt zu Beschädigungen an den Türen oder der Module. Überprüfen Sie, dass die Tragegriffe korrekt angeschraubt sind und nicht herausrutschen können.



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.

ACHTUNG

Selbst wenn Sie den Vanquish-Systemturm nur geringfügig auf dem Arbeitstisch verschieben, kann der Systemsockel beschädigt und dessen Funktionalität beeinträchtigt werden.

Bevor Sie den Systemturm verschieben, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln.

Hinweise für das Erweitern des Systemturms

Beachten Sie Folgendes:



WARNUNG—Kippgefahr beim Systemturm

Eine Konfiguration mit nur einem Turm kann die maximal erlaubte Höhe überschreiten. Wenn die Systemkonfiguration höher ist, stellt der Systemturm eine Kippgefahr dar, von der eine Verletzungsgefahr ausgehen kann.

Wenn Ihre Konfiguration höher ist, müssen Sie den Turm mit einem der folgenden Sicherheitsmaßnahmen stabilisieren:

- Verwenden Sie eines der Kits zur Stabilisierung des Turms (siehe Tabelle unten).
- Bauen Sie das System in zwei separaten Türmen auf.

Die Tabelle zeigt, welches Kit zur Stabilisierung des Turms Sie für welche Systemhöhe benötigen. Die Höhen werden von der Tischoberfläche bis zur Solvent Rack-Reling gemessen. Aufgrund von Produktions- und Installations-Toleranzen kann die tatsächliche Höhe des Turms bis zu 2 cm höher sein als die maximale Höhe. Diese Toleranz ist akzeptabel.

Turmhöhe	Erforderliches Kit zur Stabilisierung
Bis zu 100 cm	Keines Beispiel: Konfiguration mit Pumpe, Autosampler, Säulenthermostat und Variablen Wellenlängendetektor
Bis zu 123 cm	Eines der folgenden Kits: • Bench Clamp Kit • Stack Stabilizer Kit • IonBench mit Montage-Kit für den Systemturm
Bis zu 139 cm	Bench Clamp Kit
139 cm und höher	Bauen Sie das System in zwei separaten Türmen auf

5.2 Anforderungen an den Aufstellungsort

Angemessene Umgebungsbedingungen sind wichtig, um den optimalen Betrieb des Systems zu ermöglichen.

Dieser Abschnitt behandelt wichtige Anforderungen an den Aufstellungsort. Beachten Sie Folgendes:

- Stellen Sie sicher, dass sich am Aufstellungsort Steckdosen in ausreichender Anzahl für alle Geräte im System befinden.
- Betreiben Sie das System nur bei angemessenen Laborbedingungen
- Informationen zur Spezifikation finden Sie in Spezifikationen (> Seite 387).

5.2.1 Arbeitstisch

Das Vanquish-System wurde für die Aufstellung auf einem Arbeitstisch entwickelt. Wenn ein komplettes Vanquish-System und ggf. weitere Geräte auf dem Arbeitstisch aufgebaut werden soll, muss der Arbeitstisch das Gewicht aller Geräte und Komponenten tragen können, einschließlich der Lösungsmittel.

Die Abmessungen und Gewichte finden Sie in Spezifikationen (> Seite 387).

Anforderungen an den Arbeitstisch

Teil	Anforderungen
Aufstellort für den Arbeitstisch	Der Arbeitstisch muss sicher und eben auf einer vibrationsfreien Oberfläche stehen.
Arbeitstisch	 Hohe Stabilität Gute Standfestigkeit (kippsicher) Eine Höhe, die es bequem ermöglicht, den Innenraum aller Geräte im System zu erreichen. Die Tischplatte sollte trocken, sauber, sowie beständig gegen Chemikalien sein.
	 Wenn der Systemturm mit dem Bench Clamp Kit stabilisiert werden soll, muss die Arbeitsplatte des Labortisches <i>eines</i> der folgenden Anforderungen erfüllen: Stärke: Zwischen 15 und 50 mm oder
	 Material: Weich genug f ür Holzschrauben

Einzuhaltende Abstände vom System



Abbildung 24: Einzuhaltende Abstände vom System (links: Vorderansicht des Systemturms, rechts: System von oben gesehen)

Nr.	Beschreibung
1	Mindestens 10 cm
2	Mindestens 30 cm
3	Mindestens 5 cm
4	Mindestens 15 cm
5	Mindestens 35 cm

5.2.2 Hinweise zur Stromversorgung

Die Netzteile der Geräte verfügen über einen weiten Eingangsspannungsbereich und können sich automatisch an jede Netzspannung innerhalb des spezifizierten Bereichs für die Geräte 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.

Das Gerät kann durch kurze Spannungsunterbrechungen beeinträchtigt werden, wenn die Stromversorgung des Gerätes ausgeschaltet wird, um das Gerät vor einem unbestimmten Verhalten zu schützen. Bei Rückkehr der Versorgungsspannung startet die Stromversorgung das Gerät automatisch neu und eine erneute Verbindung zum Chromatographie-Datensystem ist nötig.

5.2.3 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.

5.2.4 Kondensation

ACHTUNG—Kondensation im Geräteinneren kann die Elektronik 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.

5.2.5 Betriebsbedingungen

	Stellen Sie sicher, dass der Aufstellungsort beim Betrieb des Systems die folgenden allgemeinen Umgebungs- und Betriebsbedingungen erfüllt. Angaben zur Umgebungstemperatur und Luftfeuchtigkeit finden Sie in Spezifikationen (> Seite 387).
Temperatur	
	Temperaturschwankungen können die Systemleistung beeinflussen. Vermeiden Sie signifikante Temperaturschwankungen und Luftzug. Stellen Sie das System beispielsweise nicht ins direkte Sonnenlicht, in die Nähe von Heiz- oder Kühlquellen, oder unter einen Lüftungsschacht.
Luftfeuchtigkeit	
	Die relative Luftfeuchtigkeit der Betriebsumgebung ist wichtig für die Systemleistung. Betreiben Sie das System im spezifizierten Bereich, ohne Kondensation.
	Der Betrieb des Systems in einer Umgebung mit hoher Luftfeuchtigkeit kann Kondensation und damit Schäden an den elektronischen Komponenten zur Folge haben. Bei sehr geringer Luftfeuchtigkeit kann es zur Akkumulation statischer Elektrizität und zu deren Entladung kommen, was sich negativ auf die Lebensdauer der elektronischen Bauteile auswirkt.
Belüftung	
	Der Aufstellungsort muss jederzeit gut belüftet sein, damit Gefahren für Gesundheit und Sicherheit vermieden werden können, die durch den Umgang mit gefährlichen Substanzen, flüchtigen Bestandteilen oder Gasen entstehen können.

Vibrationen und mechanische Stöße

Vibrationen und mechanische Stöße können die Systemleistung beeinflussen. Der Aufstellungsort sollte daher vibrationsfrei sein. Vermeiden Sie mechanische Stöße und stellen Sie das System nicht in der Nähe von Geräten auf, die Vibrationen verursachen.

Elektromagnetische Interferenz

- Das System wurde geprüft und erfüllt die Immunitätsanforderungen sowohl für grundlegende als auch für industrielle elektromagnetische Umgebungen gemäß EN IEC 61326-1. Starke EMV-Störquellen können die Leistung des Systems beeinträchtigen. Betreiben Sie in der Nähe des Systems keine Geräte, die starke elektrische Felder oder Magnetfelder erzeugen, sondern betreiben Sie in der Nähe des Systems nur zertifizierte Laborgeräte. Schließen Sie keine Geräte, die Stromschwankungen verursachen könnten, an dasselbe Stromnetz an, das auch das System mit Strom versorgt.
- Das System wurde geprüft und erfüllt die Emissionsgrenzwerte für Geräte der Klasse A gemäß EN IEC 61326-1.



VORSICHT

Dieses Gerät ist nicht für die Verwendung in Wohnbereichen vorgesehen und bietet in derartigen Umgebungen keinen ausreichenden Schutz vor Funkstörungen.

5.3 Aufstellen des Systems

Ein Servicetechniker von Thermo Fisher Scientific installiert das Vanquish Neo-System einschließlich aller mitgelieferten Module, Optionen und Teile, und nimmt das System in Betrieb. Der Servicetechniker prüft, dass das Vanquish Neo-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 Systems durch eine andere Person als einen Thermo Fisher Scientific-Servicetechniker erfolgt, folgen Sie den Schritten in diesem Kapitel.

Die unten aufgeführten Schritte stellen die grobe Installationsprozedur dar. Führen Sie die einzelnen Schritte in den angegebenen Kapiteln in der gezeigten Reihenfolge durch.





5.4 Öffnen und Schließen der Gerätetüren

Direkt hinter den Gerätetüren sind die für den Anwender zugänglichen Komponenten und Flussverbindungen der Geräte leicht erreichbar.

Öffnen der Gerätetüren

- 1. Öffnen Sie die linke Gerätetür.
- 2. Öffnen Sie die rechte Gerätetür.

Öffnungsmechanismus des
AutosamplersDie linke Tür des Autosamplers verfügt über einen Mechanismus, der die
Tür automatisch öffnet, wenn der Autosampler ausgeschaltet ist, um
eine ordnungsgemäße Belüftung des Probenraums zu ermöglichen. Bei
einem Stromausfall, oder wenn das Netzkabel gezogen wird, während
der Autosampler einschaltet ist, öffnet der Mechanismus die linke
Gerätetür automatisch.





Schließen der Gerätetüren

- 1. Schließen Sie die rechte Gerätetür.
- 2. Schließen Sie die linke Gerätetür.

Schließmechanismus des Wenn der Autosampler eingeschaltet ist, kann die linke Gerätetür Autosamplers geschlossen werden.

5.5 Installation des Säulenthermostaten

5.5.1 Einbau des Säulenthermostaten in den Systemturm

Der Säulenthermostat ist für das Vanquish-System optional. Die Module des Systems werden in einem Systemturm aufgebaut, wobei die genaue Anordnung von der Systemkonfiguration abhängt.

Erforderliche Werkzeuge

- Schraubendreher, Torx T10
- Schraubendreher, Torx T20

Vorbereitungen

Wenn das System bereits in Betrieb ist:

- 1. Schalten Sie den Pumpenfluss aus.
- 2. Schalten Sie das System aus.
- 3. Trennen Sie die Schläuche von den Behältern.
- 4. Entfernen Sie die Behälter aus dem Solvent Rack.

Gehen Sie wie folgt vor

- 1. Lösen Sie mit einem Schraubendreher (Torx T10) die Schrauben, die die Blechklammer am Autosampler befestigen.
- 2. Entfernen Sie die Blechklammer.

TIPP Bewahren Sie die Blechklammer und die Schrauben auf. Sie benötigen diese gegebenenfalls bei einem künftigen Transport.



Abbildung 26: Entfernen der Blechklammer (hier: Autosampler)

3. Heben Sie das Solvent Rack vorne an und ziehen Sie das Solvent Rack nach vorne heraus.

- 4. Heben Sie den Säulenthermostaten mit Hilfe der Tragegriffe an. Setzen Sie den Säulenthermostaten ca. 5 cm vor dem Ende der Schiene auf dem Autosampler auf. Schieben Sie den Säulenthermostaten nach hinten, bis er einrastet.
- 5. Lösen Sie die Schrauben an den Tragegriffen mit einem Schraubendreher (Torx T20), bis Sie den Tragegriff in der Schiene bewegen können. Entfernen Sie die Schrauben nicht vollständig von den Tragegriffen.



Abbildung 27: Tragegriffe

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

6. Ziehen Sie die Tragegriffe zur Rückseite des Säulenthermostaten hin aus den Schienen heraus.



Abbildung 28: Herausziehen des Tragegriffs aus der linken Schiene

- 7. Setzen Sie das Solvent Rack ca. 5 cm vor dem Ende der Schiene auf dem Säulenthermostaten auf. Schieben Sie das Solvent Rack nach hinten, bis es einrastet.
- 8. Verbinden Sie die Schläuche mit den Behältern.
- 9. Schließen Sie den Drainage-Ablauf des Säulenthermostaten an die Systemdrainage an.
- 10. Schließen Sie das Netzkabel, die USB- und System-Interlink-Kabel an den Säulenthermostaten an.

Sehen Sie dazu auch

- Anschließen der Netzkabel (> Seite 115)
- Anschließen der USB-Kabel (> Seite 109)

- Anschließen der System-Interlink-Kabel (> Seite 113)
- Kapillar- und Schlauchführung durch das System (> Seite 119)

5.5.2 Installieren des Vanquish-Schaltventils

Sie können Vanquish-Schaltventile an der rechten oder linken Position installieren, wenn kein VICI-Ventilantrieb installiert ist.

Erforderliche Teile

Vanquish-Schaltventil

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Vorbereitungen

- 1. Schalten Sie den Säulenthermostaten aus. Wenn der Säulenthermostat eingeschaltet ist, kann das Schaltventil während der Installation beschädigt werden.
- 2. Entfernen Sie die Verschlussfittinge noch nicht von den Ventilports.

Gehen Sie wie folgt vor

- Um die Abdeckung zu entfernen, schrauben Sie einen Verschlussfitting von den Ventil-Ports in die Öffnung in der Mitte der Abdeckung.
- 2. Lösen Sie die 4 Schrauben an der Abdeckung mit dem Schraubendreher.
- 3. Ziehen Sie die Abdeckung mit dem Verschlussfitting heraus.



Abbildung 29: Entfernen der Abdeckung



4. Entfernen Sie das Klebeband, das den Kabelstecker abdichtet.

Abbildung 30: Entfernen des Klebebands

5. Ziehen Sie das Kabel aus dem Gehäuse heraus und schließen Sie das Kabel an den Stecker auf der Rückseite des Schaltventils an.



Abbildung 31: Anschließen des Kabels an den Stecker

- 6. Schieben Sie das Kabel zurück ins Gehäuse.
- 7. Schieben Sie das Schaltventil vorsichtig ins Gehäuse hinein.



Abbildung 32: Installieren des Vanquish-Schaltventils

8. Prüfen Sie die Ausrichtung des Schaltventils: Port 3 muss nach oben zeigen.

9. Ziehen Sie die 4 Schrauben am Schaltventil fest.

Abbildung 33: 4 Schrauben am Schaltventil

- 10. Entfernen Sie die Verschlussfittinge von den Ventilports.
- 11. Schalten Sie den Säulenthermostaten ein.
- 12. Führen Sie das erforderliche Skript aus, um Ihren Workflow zu ändern, und folgen Sie den Schritten im Skript, um die Schläuche zu verbinden.
- 13. Prüfen Sie die Ventile visuell auf Undichtigkeit.
- Führen Sie das Skript aus: D02 Systemdichtigkeit testen (was getestet werden soll = System). Dieses Skript testet auch die Säulenschaltventile auf Undichtigkeit.

5.5.3 Installieren des VICI-Schaltventils

Sie können VICI-Schaltventile nur dort installieren, wo ein VICI-Ventilantrieb installiert ist. Um den VICI-Ventilantrieb zu installieren, wenden Sie sich an den Kundendienst von Thermo Fisher Scientific. Eine Übersicht des VICI-Schaltventils finden Sie unter Übersicht der Säulenschaltventile () Seite 54).

Voraussetzung

Der VICI-Ventilantrieb und der Adapterrahmen sind installiert (durch Servicetechniker erfolgt).

Erforderliche Teile

VICI-Schaltventil

Erforderliche Werkzeuge

• Schraubendreher, Torx T10

Gehen Sie wie folgt vor

- Um die Abdeckung zu entfernen, schrauben Sie einen Verschlussfitting von den Ventil-Ports in die Öffnung in der Mitte der Abdeckung.
- 2. Lösen Sie die 4 Schrauben an der Abdeckung mit dem Schraubendreher.
- 3. Ziehen Sie die Abdeckung mit dem Verschlussfitting heraus.



Abbildung 34: Entfernen der Abdeckung

- 4. Entfernen Sie die Verschlussfittinge noch nicht von den Ventilports.
- 5. Richten Sie die beiden Stifte auf der Rückseite des Ventilklemmrings an den entsprechenden Öffnungen im Ventilflansch aus.



Abbildung 35: Ausrichten des Ventilkopfes am Flansch

- 6. Drücken Sie den Ventilkopf auf den Ventilflansch.
- Um die Achse mit dem Ventilkopf auszurichten, halten Sie je nach Ventilposition die Taste VALVE LEFT oder VALVE RIGHT auf der Tastatur gedrückt. Die Funktion "Catch Valve" wird aktiviert und Sie können die Taste loslassen.
- 8. Wenn die Achse ausgerichtet ist, rastet der Ventilkopf in den Ventilflansch ein.

- 9. Entfernen Sie die Verschlussfittinge von den Ventilports.
- 10. Stellen Sie die Verbindungen je nach Erfordernis für Ihre Anwendung her.
- 11. Konfigurieren Sie das VICI-Schaltventil im Chromatographie-Datensystem.
- 12. Prüfen Sie die Ventile visuell auf Undichtigkeit.
- Führen Sie das Skript aus: D02 Systemdichtigkeit testen (was getestet werden soll = Sampler). Dieses Skript testet auch die Säulenschaltventile auf Undichtigkeit.

5.5.4 Installieren des passiven Vorheizers

Der passive Vorheizer ist optional. Er kann verwendet werden, um den Eluenten an die Säulentemperatur anzupassen, bevor der Eluent die Säule erreicht.

Der passive Vorheizer ist für Flussraten > 50 μ L/min, insbesondere für Säulen mit einem Innendurchmesser von 1 mm, in einem Flussbereich von 50-100 μ L/min gedacht.

Erforderliche Teile

- Passiver Vorheizer
- 2 Schrauben

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Gehen Sie wie folgt vor

 Positionieren Sie den passiven Vorheizer an der gewünschten Stelle im Säulenthermostaten.
 Im Säulenraum befinden sich zwei verschiedene Einbaupositionen.

TIPP

Der passive Vorheizer kann an jeder der zwei unten dargestellten Stellen positioniert werden. Die Ausrichtung des passiven Vorheizers kann je nach Erfordernis erfolgen: die Auslasskapillare kann sowohl nach unten als auch nach oben zeigen, je nach Position des Säuleneingangs.



Abbildung 36: Einbaupositionen für den passiven Vorheizer

- 2. Richten Sie die Schraublöcher des passiven Vorheizers und die Schraublöcher an der jeweiligen Einbauposition aneinander aus.
- 3. Ziehen Sie die zwei Schrauben und Unterlegscheiben mit einem Schraubendreher fest, um den passiven Vorheizer im Säulenraum zu befestigen.
- 4. Schließen Sie die Auslasskapillare des passiven Vorheizers an den Säuleneingang an.
- Schließen Sie die Einlasskapillare des passiven Vorheizers an den Autosampler oder ein Säulenschaltventil an. Wenn Sie den passiven Vorheizer an den Autosampler anschließen, führen Sie die Kapillare durch die untere Kapillarführung.



Abbildung 37: Kapillarführung für das Anschließen an den Autosampler

ΤΙΡΡ

Die Temperatur des passiven Vorheizers wird nicht geregelt. Daher ist es nicht erforderlich den passiven Vorheizer im Chromatographie-Datensystem einzurichten.

5.5.5 Installieren von Säulen im Säulenthermostaten

Im Säulenraum sind acht Säulenhaltespangen installiert. Befestigen Sie die Säulen an den Säulenhaltespangen mit Hilfe der Befestigungsclips aus dem Zubehörkit. Sie können die Befestigungsclips in jeder beliebigen Position an den Säulenhaltespangen anbringen.



Abbildung 38: Mögliche Positionen der Säulenbefestigungsclips

Erforderliche Teile

- Säule, die für Ihre Anwendung erforderlich ist
- Säulenbefestigungsclips
- Selbstklebende Schaumstoffpads

Gehen Sie wie folgt vor

1. Drücken Sie den Ring in Pfeilrichtung leicht zusammen, um ihn vom Unterteil des Säulenbefestigungsclips abzuziehen.



Abbildung 39: Säulenbefestigungsclips

Nr.	Beschreibung
1	Ring
2	Schaumstoffpad
3	Unterteil

- 2. Im Lieferumfang der Befestigungsclips sind selbstklebende Schaumstoffpads enthalten. Kleben Sie zum Schutz der Säule auf der Innenseite jedes Unterteils eines dieser Pads auf.
- Setzen Sie das Unterteil des S\u00e4ulenbefestigungsclips an der gew\u00fcnschten Stelle auf die Haltespange auf und drehen Sie es um 90° im Uhrzeigersinn.
- 4. Stellen Sie die Kapillarverbindungen an Säuleneingang und Säulenausgang her.
- 5. Legen Sie die Säule in das Unterteil des Befestigungsclips und schieben Sie den Ring auf.



Abbildung 40: Aufschieben des Rings

6. Führen Sie die Kapillaren durch eine der in der Abbildung gezeigten Kapillarführungen.



Abbildung 41: Kapillarführungen

- Um die S\u00e4uleneigenschaften einzustellen, f\u00fchren Sie die Skripte aus: A03 - Trenns\u00e4ulentyp festlegen und A04 -Trenns\u00e4ulenspezifikationen festlegen (und A05 – Trap-S\u00e4ulenspezifikationen festlegen nach Bedarf).
- 8. Schalten Sie den Pumpenfluss ein.

5.6 Entfernen der Transportsicherungen des Autosamplers

Erforderliche Werkzeuge

Schlitzschraubendreher

Gehen Sie wie folgt vor

 Entfernen Sie sonstige Transportsicherungen oder Schutzabdeckungen vom Autosampler wenn vorhanden. Transportsicherungen sind entsprechend zum Entfernen markiert.



2. Öffnen Sie die Autosamplertüren.

Abbildung 42: Öffnen der Autosamplertüren

 Entfernen Sie die Isolierabdeckung (siehe Entfernen der Isolierabdeckung (▶ Seite 98)).
 Bewahren Sie die Abdeckung für den Betrieb auf.

ACHTUNG

Wenn die Transportsicherungsschraube nicht korrekt gelöst wurde, kann der Autosampler beschädigt werden. Vergewissern Sie sich, dass die Transportsicherungsschraube vollständig gelöst ist, bevor Sie den Autosampler einschalten.

 Lösen Sie die Transportsicherungsschraube oberhalb der Nadeleinheit: Drehen Sie die unverlierbare Schraube gegen den Uhrzeigersinn mit Ihren Fingern, bis sie locker in ihrer Feder sitzt. Die Verwendung eines Schlitzschraubendrehers erleichtert gegebenenfalls das Drehen der Schraube.





- Montieren Sie die Isolierabdeckung (siehe Einbauen der Isolierabdeckung (▶ Seite 99)).
- 6. Entfernen Sie den Schnappniet von der Drehteller-Transportsicherung und entnehmen Sie die Drehteller-Transportsicherung.

5.6.1 Entfernen der Isolierabdeckung

Gehen Sie wie folgt vor

- 1. Nehmen Sie die Abdeckung am unteren Ende und ziehen Sie sie leicht, um die Isolierabdeckung von der Halteklammer zu lösen.
- 2. Ziehen Sie die Abdeckung vorsichtig, um sie aus dem oberen Gehäuserahmen herauszuziehen.
- 3. Ziehen Sie die Abdeckung von der Halteklammer weg.



Abbildung 44: Entfernen der Isolierabdeckung

5.6.2 Einbauen der Isolierabdeckung

Erforderliche Teile

Isolierabdeckung

Gehen Sie wie folgt vor

1. Legen Sie die Probenschleife richtig in den dafür vorgesehenen Haken an der Halteklammer der Isolierabdeckung ein.



Abbildung 45: Isolierabdeckung mit Halteklammer (links); Halteklammer mit Haken für die Probenschleife (rechts)

- 2. Richten Sie die Abdeckung an der Nadeleinheit und an der Probenschleife an der inneren Frontblende aus.
- 3. Um die Abdeckung einzubauen, drücken Sie die Nut in der Abdeckung vorsichtig in die Halteklammer. Achten Sie darauf, dass die Probenschleife nicht durch die Abdeckung gequetscht wird.

5.7 Montieren des Displays

Sie können das Display an der linken Seite von jedem Systemturmmodul montieren (wenn das Modul den Anforderungen unter *Voraussetzung* entspricht). Um das Display zu drehen, halten Sie auf der linken Seite einen Abstand von 10 cm ein.

Voraussetzung

Sie können das Display nur an Module montieren, die eine kantige Aussparung besitzen. Das Display kann nicht an Gehäusen mit einer Uförmigen Aussparung montiert werden.



Abbildung 46: Gehäuse mit Aussparungen (links: eckig, rechts: U-förmig)

Vorbereitungen

Packen Sie das Display aus.

TIPP Bewahren Sie die Originalverpackung und alle Verpackungsmaterialien auf. Sie benötigen diese Dinge, wenn Sie das Display an einen anderen Ort transportieren möchten.

Erforderliche Teile

- Display-Baugruppe mit:
 - ♦ 3 Torx[™]-Schrauben
 - Nutenstein
 - Gleitnutenstein



Abbildung 47: Gleitnutenstein und Nutenstein

Nr.	Beschreibung
1	Gleitnutenstein
2	Nutenstein

- Display-Sperre
- Kabelabdeckung

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Gehen Sie wie folgt vor

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

1. Prüfen Sie, ob das Display-Kabel wie unten gezeigt verläuft.



Abbildung 48: Korrekter Verlauf des Display-Kabels

Nr.	Beschreibung
1	Schwenkarm
2	Display-Kabel

2. Wenn das Display-Kabel nicht wie in der Abbildung verläuft, drehen Sie das Display um den Schwenkarm im bzw. gegen den Uhrzeigersinn bis das Display-Kabel so verläuft wie oben gezeigt.



Abbildung 49: Drehen des Displays um den Schwenkarm

3. Lockern Sie die oberen 2 Schrauben und drehen Sie die untere Schraube am Display heraus, um den Nutenstein zu entfernen.



Abbildung 50: Betroffene Schrauben am Display

4. Schieben Sie den Gleitnutenstein des Displays von hinten in die linke Schiene des Moduls und ziehen Sie das Display bis zur Sperre nach vorne.



Abbildung 51: Ziehen des Displays

Nr.	Beschreibung
1	Display
2	Gleitnutenstein
3	Schiene



5. Ziehen Sie die 2 Schrauben in den oberen zwei Bohrungen etwas an.

Abbildung 52: Befestigen des Displays

6. Schauen Sie nach der Aussparung auf der Innenseite der linken Gehäusewand.



Abbildung 53: Aussparung

7. Halten Sie den Nutenstein wie gezeigt auf der Höhe der Aussparung. Halten Sie den Nutenstein, so dass die Bohrungen im Nutenstein und im Gehäuse fluchten.





 Halten Sie den Nutenstein mit der rechten Hand und befestigen Sie ihn leicht mit der linken Hand von außen mit einer Schraube.
 Benutzen Sie dabei eine der Bohrungen, die in der Abbildung unten mit einem Kreis gekennzeichnet sind.



Abbildung 55: Befestigen des Nutensteins

- 9. Prüfen Sie, ob sich das Display in horizontaler Position befindet. Wenn nicht, korrigieren Sie die Ausrichtung des Displays wie folgt:
 - a) Bewegen Sie das Display bis es waagrecht ist.
 - b) Ziehen Sie die 3 Schrauben fest an.
- 10. Wenn Sie das Display an das oberste Systemmodul montiert haben, klemmen Sie die Display-Sperre wie gezeigt an das Solvent Rack.



Abbildung 56: Display-Sperre

5.8 Anschließen der Kabel

5.8.1 Übersicht der Anschlüsse an den Systemmodulen

Folgende Anschlüsse stehen an den Systemmodulen für Netzanschluss und Signalanschlüsse zur Verfügung:



Abbildung 57: Anschlüsse

Nr.	Beschreibung
1	System Interlink-Port Für das Ein- und Ausschalten des Systems über den Vanquish-Systemsockel und Kommunikation zwischen den Geräten Siehe Anschließen der System-Interlink-Kabel (▶ Seite 113).
2	Digital I/O-Ports (Dig I/O) Für den Austausch digitaler Signale mit externen Geräten. Verwenden Sie Dig I/O 2 am Autosampler, um ein Massenspektrometer anzuschließen. Siehe Anschließen eines Massenspektrometers (▶ Seite 117).
3	USB-Hub (Stecker Typ "A") Für den Anschluss anderer Module des Vanquish-Systems
4	USB-Port (Stecker Typ "B") Für den Anschluss an andere Module im Vanquish-System oder an den Rechner, auf dem das Chromatographie-Datensystem installiert ist Siehe Anschließen der USB-Kabel () Seite 109).
5	Typenschild (Beispiel) mit Angabe von Seriennummer, Teilenummer, Modulnamen, Revisionsnummer, Netzspannung und Nennleistung der Sicherungen
6	Hauptnetzschalter (Ein/Aus)
7	Sicherungshalter
8	Netzbuchse

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.

5.8.2 Übersicht der Anschlüsse am System Controller

Folgende Anschlüsse befinden sich auf der kurzen Seite des System Controllers:



Abbildung 58: Anschlüsse auf der kurzen Seite

Nr.	Beschreibung
1	Typenschild
	Gibt Auskunft über Seriennummer und Netzspannung.
2	Vanquish System Controller (VSC)-Port
	Für den Anschluss an den Systemsockel.
3	System Interlink-Port
	Für den Anschluss an den Systemsockel und an ein Modul im HPLC-System
4	USB-Port (Universal Serial Bus) (Stecker Typ "A")
	Deaktiviert - Nicht verwenden
5	Local Area Network (LAN)-Port
	Für den Anschluss an ein firmeninternes Netzwerk.
6	Hauptnetzschalter (Ein/Aus)
7	Netzbuchse
Die folgenden Anschlüsse und Steckplätze befinden sich auf der langen Seite des System Controllers:



Abbildung 59: Anschlüsse auf der langen Seite

Nr.	Beschreibung
1	Netzbuchse für das Display
2	Datensignal-Port für das Display
3	System Interlink-Ports Deaktiviert
4	USB-Port (Stecker Typ "A") Deaktiviert - Nicht verwenden
5	Schild mit Angabe von Teilenummer, Modulnamen und Revisionsnummer.
6	Storage Device (SD)-Kartensteckplatz

5.8.3 Anschließen der USB-Kabel

Für alle USB-Verbindungen wird ein Standard-USB-Kabel (Typ A auf Typ B) benötigt.



Abbildung 60: USB-Kabel

Der flache, rechteckige Stecker ist der Stecker Typ "A". Der kleinere, sechseckige Stecker ist der Stecker Typ "B".

TIPP

- Ein USB-Kabel darf laut Norm maximal 5 m lang sein. Die Länge der USB-Verbindung zum Rechner oder nächsten USB-Hub darf 5 m nicht überschreiten.
- Vergewissern Sie sich nach dem Anschluss der USB-Kabel, jedoch bevor Sie ein Gerät zum ersten Mal einschalten, dass die Chromatographie-Software auf dem Rechner installiert ist. Die benötigten USB-Treiber werden automatisch geladen und das Windows[™]-Betriebssystem kann das Gerät erkennen, wenn dieses eingeschaltet ist.



Abbildung 61: USB-Verbindungen (hier: mit installiertem Säulenthermostat)

Nr.	Beschreibung
-	USB-Anschluss:
1	Von der Pumpe zum Rechner
2	Vom Autosampler zur Pumpe
3	Vom Säulenthermostat (falls installiert) zur Pumpe

Gehen Sie wie folgt vor

- Schließen Sie an Autosampler, Pumpe und Säulenthermostat (falls installiert) das USB-Kabel mit dem Stecker Typ "B" an den Port USB an.
- 2. Schließen Sie das andere Ende des USB-Kabels von der Pumpe (Typ "A") an den Hub **USB** (2.0 oder höher) am Rechner an.
- 3. Schließen Sie das andere Ende des USB-Kabels von den anderen Modulen (Typ "A") an den Hub **USB** an der Pumpe an.

Eine Steuerung der Module, die über einen Hub **USB** an einem anderen Modul angeschlossen sind, ist nur möglich, wenn das Modul mit dem Hub eingeschaltet ist.

ACHTUNG

- Verwenden Sie keine beschädigten Kommunikationskabel. Tauschen Sie das Kabel aus, wenn Sie den Verdacht haben, das Kabel sei beschädigt.
- Verwenden Sie nur die von Thermo Fisher Scientific f
 ür den Anschluss des Systems angebotenen Anschlusskabel, um eine einwandfreie Funktion der Verbindung zu gew
 ährleisten.

5.8.4 Anschließen der Display-Kabel

Erforderliche Werkzeuge

Schere, um die Kabelabdeckung zurecht zu schneiden

Gehen Sie wie folgt vor

- 1. Verbinden Sie das Display-Netzkabel mit dem **Display DC OUT 24V**-Port am System Controller.
- Befestigen Sie das Display-Datensignalkabel fingerfest am Display Data-Port des System Controllers.



3. Verlegen Sie das Display-Netzkabel und das Datensignalkabel in der Schiene.

Abbildung 62: Verlegen des Display-Netzkabels

- Kürzen Sie die Kabelabdeckung mit der Schere auf eine Länge von 34 cm.
- 5. Drücken Sie die Kabelabdeckung fest in die Schiene.



Abbildung 63: Befestigen der Kabelabdeckung



6. Fixieren Sie den senkrechten Teil des Kabelstrangs in den Kabelhalterungen.

Abbildung 64: Fixieren der Kabel mit den Kabelhalterungen

- 7. Schieben Sie den Systemturm zurück in Betriebsposition.
- 8. Sperren Sie den Systemsockel mit den Systemsockelschlüsseln (siehe Sperren des Systemsockels (▶ Seite 118)).

5.8.5 Anschließen der System-Interlink-Kabel

Für die System-Interlink-Anschlüsse ist ein spezielles System-Interlink-Kabel erforderlich (Cat. 6-Kabel).



Abbildung 65: System-Interlink-Kabel

Die System-Interlink-Verbindungen von einem Systemmodul zum anderen sind bis auf die Verbindungen zum optionalen Säulenthermostat vorinstalliert.



Abbildung 66: System-Interlink-Anschlüsse

Nr.	Beschreibung
-	System-Interlink-Verbindung:
1	Vom Systemsockel zum System-Controller
2	Vom System-Controller zur Pumpe
3	Von der Pumpe zum Autosampler
4	Vom Autosampler zum Säulenthermostat (falls installiert)

Wenn ein Säulenthermostat installiert ist (optional)

- 1. Entfernen Sie die Kappe, die den rechten Port **System Interlink** am Autosampler abdeckt.
- Schließen Sie das eine Ende des System-Interlink-Kabels an den Port System Interlink am Säulenthermostaten an, und verbinden Sie das andere Ende mit dem rechten Port System Interlink am Autosampler (wie in der Abbildung gezeigt).

5.8.6 Anschließen eines LAN-Kabels

Erforderliche Teile

- LAN-Kabel
- Netzwerkkarte

Gehen Sie wie folgt vor

ACHTUNG—Schäden am Vanquish-Gerät

Die **System Interlink**- und die **LAN**-Ports besitzen denselben Steckertyp. Wenn Sie zum Beispiel ein LAN-Kabel an einen **System Interlink**-Port eines Vanquish-Moduls anschließen, kann dies das Vanquish-Modul und den System Controller beschädigen.

Verbinden Sie nur Ports desselben Schnittstellentyps.

- 1. Schließen Sie ein LAN-Kabel an den LAN-Port des System Controllers an.
- Schließen Sie die andere Seite des LAN-Kabels an den LAN-Port der Netzwerkkarte im Datensystemrechner, auf dem die Chromatographie-Software installiert ist.

5.8.7 Anschließen der Netzkabel

In diesem Abschnitt finden Sie Einzelheiten darüber, wie Sie die Module an die Stromversorgung anschließen. Schalten Sie System und Module noch nicht ein.

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.

Gehen Sie wie folgt vor

- 1. Vergewissern Sie sich, dass alle Geräte über den Netzschalter ausgeschaltet sind.
- 2. Verbinden Sie das Netzkabel an allen Geräten mit der Netzbuchse.
- 3. Verbinden Sie das andere Ende der Netzkabel mit einer geeigneten Stromquelle.
- 4. Stecken Sie die Signal- und Netzkabel in die Kabelhalterungen (siehe Verlegen der Signalkabel in den Kabelhalterungen (▶ Seite 116)).
- 5. Wenn nötig, bringen Sie das System an seine endgültige Position.
- 6. Sichern Sie den Systemsockel gegen Verschieben (siehe Sperren des Systemsockels (▶ Seite 118)).

5.8.8 Verlegen der Signalkabel in den Kabelhalterungen

Mit den Kabelhalterungen an der Systemrückseite können Sie Signalkabel ordentlich am System entlangführen. Die Pumpe und der optionale Säulenthermostat besitzen jeweils eine Kabelhalterung auf jeder Seite und der Autosampler zwei Kabelhalterungen auf jeder Seite. Das Bild zeigt, wo sich die Kabelhalterungen befinden.



Abbildung 67: Kabelhalterungen am System

5.9 Anschließen eines Massenspektrometers

Erforderliche Teile

Digitales I/O-Signalkabel



Abbildung 68: Digitales I/O-Signalkabel

Pin	Beschreibung
1	6-poliger Stecker, für die Verbindung zum Dig I/O-Port am System
2	6 Adern, für die Verbindung mit einem externen Gerät, zum Beispiel, einem Massenspektrometer

Gehen Sie wie folgt vor

- 1. Stecken Sie am Autosampler den 6-poligen Stecker des Kabels in den Port mit der Kennzeichnung **Dig I/O 2**.
- Verbinden Sie die Adern f
 ür die Ausgangssignale (weiße Ader und braune Ader) mit den geeigneten Anschl
 üssen am Massenspektrometer.
 Damit wird das Inject-Response-Signal automatisch
 übertragen.

5.10 Sperren des Systemsockels

Nach dem Auspacken des Systems wurde es in den Verschiebemodus versetzt. Deswegen stecken die Systemsockelschlüssel bereits im Systemsockel.

- 1. Schieben Sie den Systemturm in Betriebsposition.
- 2. Sperren Sie den Systemsockel, indem Sie die Schlüssel wie unten gezeigt drehen.



Abbildung 69: Sperren des Systemsockels (Ansicht von oben)

3. Ziehen Sie die Schlüssel ab und bewahren Sie sie an einem sicheren Platz auf, zum Beispiel in der Schublade des Systemsockels.

5.11 Anschließen der Flussverbindungen

Bei der Auslieferung des Systems sind die meisten Flussverbindungen für den Workflow "Direkte Injektion" vorinstalliert. Folgen Sie den Anweisungen in diesem Abschnitt, um die Installation aller Flussverbindungen abzuschließen.

TIPP Um auf einen anderen Workflow zu wechseln, schließen Sie die Installation wie in dieser Anleitung beschrieben ab und führen Sie das Skript aus: **A06 - Fluidiksystem/Workflow ändern**. Weitere Informationen zu anderen Workflows finden Sie im *Vanquish Neo System User Guide*.

5.11.1 Kapillar- und Schlauchführung durch das System

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

Kanal für die Durchführung der Schläuche, einschließlich Schlauchführungen

Um Schläuche und Verbindungen vom obersten Modul zum untersten Modul durch den Vanquish Neo-Systemturm zu führen, sind die stapelfähigen Module innen rechts mit einem Schlauchkanal ausgestattet. Der Schlauchkanal hat vier Schlauchführungen.

Drücken Sie die Schläuche oder Verbindungen in die Schlauchführungen.



Abbildung 70: Schlauchkanal mit Schlauchführungen (links: Ansicht von innen, rechts: von oben)

Schlauchklammern

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 71: Schlauchklammer (links); Schlauchklammer installiert (rechts)

Führungslöcher und Kapillarclips

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

5.11.2 Anschließen von Fittingen, Kapillaren und Schläuchen

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

5.11.2.1 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.

5.11.2.2 Anschließen von NanoViper-Kapillaren

Dieser Abschnitt beschreibt den Anschluss von nanoViper[™]-Kapillaren (im Weiteren als Viper-Kapillaren[™] bezeichnet).





Die Viper-Fittinge werden ohne Rändelschraube installiert. Bei Bedarf können Sie die Rändelschrauben verwenden, die im Zubehörkit mitgeliefert werden.

ACHTUNG

- Lösen und ziehen Sie die Viper-Kapillaren nur mit den Fingern und dem speziellen Drehmoment-Werkzeug fest. Verwenden Sie ausschließlich die für diesen Zweck von Thermo Fisher Scientific zur Verfügung gestellten Werkzeuge. Verwenden Sie das Werkzeug nur für die Verbindung, für die es vorgesehen ist.
- Um Schäden an den Kapillaren und Anschlüssen zu vermeiden, lösen und ziehen Sie die Viper-Kapillaren nur fest, wenn der Systemdruck gleich Null ist.

Erforderliche Werkzeuge

Drehmoment-Schraubendreher mit:

- Basis
- Verlängerung
- Viper-Steckaufsatz

Gehen Sie wie folgt vor

ACHTUNG—Beschädigung am nanoViper-Fitting

Fassen Sie den Drehmoment-Schraubendreher an der Basis und halten Sie ihn gerade. Durch ein Neigen des Drehmoment-Schraubendrehers wird zu viel Kraft auf das nanoViper-Fitting ausgeübt und das Fitting oder der Drehmoment-Schraubendreher können beschädigt werden.

- 1. Führen Sie die Viper-Kapillare in den Anschlussport ein.
- 2. Ziehen Sie die Verbindung mit den Fingern fest.
- Ziehen Sie die Verbindung mit dem Drehmoment-Schraubendreher weiter fest. Verwenden Sie die Verlängerung, um schwer erreichbare Kapillaren festzuziehen.



Abbildung 73: Drehmoment-Schraubendreher

Nr.	Beschreibung
1	Viper-Steckaufsatz
2	Verlängerung
3	Basis

- 4. Ziehen Sie die Verbindung fest, bis die Drehmomentbegrenzung auslöst (erkennbar am hörbaren Klicken).
- Prüfen Sie die Verbindung auf Undichtigkeit. Wenn eine Undichtigkeit vorhanden ist, beheben Sie die Undichtigkeit.

Beheben von Undichtigkeiten

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

5.11.3 Anschließen der Waste-Leitungen

5.11.3.1 Übersicht der Waste-Leitungen



WARNUNG—Austritt gefährlicher Substanzen

Gefährliche Substanzen können aus Waste-Leitungen und Abfallbehältern austreten, wenn die Verbindungen und die Abfallbehälter nicht korrekt installiert sind. Der Kontakt mit diesen Substanzen kann Gesundheits- und Sicherheitsrisiken darstellen.

Überprüfen Sie die ordnungsgemäße Installation der Verbindungen und Behälter, indem Sie das Drainagesystem testen (siehe Testen des Drainagesystems (▶ Seite 127)).

Tritt im System eine Undichtigkeit auf, fließt die Flüssigkeit über den Drainage-Ablauf am Vanquish-Systemsockel in den Abfall.

Ableiten von Flüssigkeiten in den Abfall

Diese Flüssigkeiten werden durch den Drainage-Ablauf rechts unten an der Leckage-Wanne in die Systemdrainage des Vanquish-Systems geleitet:

- Flüssigkeiten, die sich in den Leckage-Wannen sammeln
- Abfallflüssigkeiten aus den Nadel-Waschsystemen und der Drainagepumpe des Autosamplers
- Abfallflüssigkeit vom Flowmeter



Abbildung 74: Übersicht der Waste-Leitungen

Nr.	Beschreibung
1	Ablaufschlauch des Nadel-Waschsystems
2	Ablaufschlauch der Drainagepumpe
3	Flowmeter-Abfallleitung (Purge Out)

Für das Ableiten von Flüssigkeiten in den Abfall, siehe Anschließen des Ablaufschlauchs am System-Drainage-Ablauf (▶ Seite 125).

Um das Drainage-System zu testen, wenn die Waste-Leitung angeschlossen ist, siehe Testen des Drainagesystems (> Seite 127).

TIPP Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt und stellen Sie sicher, dass der Schlauch frei von Kerben und Graten ist.

5.11.3.2 Anschließen des Ablaufschlauchs am System-Drainage-Ablauf

Erforderliche Teile

- Ablaufschlauch aus dem Zubehör
- L-Stück aus dem Zubehör
- Geeigneter Abfallbehälter

Gehen Sie wie folgt vor

1. Schieben Sie ein L-Stück bis zum Anschlag in den System-Drainage-Ablauf; die Verbindungen sind dann selbstdichtend.



Abbildung 75: Verbundener Ablaufschlauch mit L-Stück

Nr.	Beschreibung
1	Trichter für Systemdrainage
2	Detektor-Waste-Auslass
2a	Detektor-Waste-Leitung
3	L-Stück
3a	System-Ablaufschlauch

- 2. Schieben Sie den Ablaufschlauch bis zum Anschlag in das L-Stück ein; die Verbindungen sind dann selbstdichtend.
- 3. Führen Sie das andere Ende des Ablaufschlauchs in einen Abfallbehälter.

4. *Gilt nur, wenn Sie den Ablaufschlauch für die Systemdrainage nach hinten führen:* Setzen Sie den Ablaufschlauch für die Systemdrainage in die Klemmen ein und stellen Sie sicher, dass der Ablaufschlauch keinen Siphon bildet.



Abbildung 76: Ablaufschlauch in Klemmen einsetzen

5. Prüfen Sie, ob die Flüssigkeit ungehindert abfließt (siehe nächster Abschnitt).

Prüfen Sie, ob die Flüssigkeit ungehindert abfließt:

- Stellen Sie den Abfallbehälter unterhalb des Systemturmniveaus auf. Damit keine Lösungsmitteldämpfe zurück ins System gelangen können, vergewissern Sie sich, dass der Abstand ausreichend groß ist.
- 2. Stellen Sie sicher, dass der Ablaufschlauch an keiner Stelle geknickt oder eingeklemmt ist.
- 3. Der Ablaufschlauch muss sich vollständig unterhalb des System-Drainage-Ablaufs befinden, damit die Flüssigkeit nicht zurück ins System fließen kann. Vermeiden Sie die Bildung von Siphons.
- 4. Damit die Abfallflüssigkeit korrekt abfließt und eine Ansammlung im Systemsockel verhindert wird, darf das Ende des Ablaufschlauchs nie in die Abfallflüssigkeit eintauchen (siehe Abbildung unten):
 - a) Kürzen Sie den Ablaufschlauch entsprechend, falls erforderlich.
 - b) Beobachten Sie den Füllstand im Abfallbehälter und entleeren Sie den Abfallbehälter, wenn erforderlich.



Abbildung 77: Schlauchende im Abfallbehälter

5.11.3.3 Testen des Drainagesystems

Prüfen Sie, ob im Drainagesystem die Flüssigkeit ungehindert abfließt:

- 1. Gießen Sie Wasser oder Isopropanol in den Ablauf des Solvent-Racks.
- 2. Prüfen Sie, ob die Flüssigkeit am System-Drainage-Ablauf austritt und in den Abfallbehälter abfließt.
- 3. Tritt an einer Stelle Flüssigkeit aus, beheben Sie die Ursache für die Undichtigkeit.

5.12 Füllen der Lösungsmittelbehälter

Füllen Sie vor dem Betrieb des Systems die Lösungsmittelbehälter.

TIPP Um die Bildung von kleinen Luftblasen in den Schläuchen zu vermeiden, platzieren Sie die Behälter im Solvent Rack oder oberhalb des Systemturms.

Vorbereitungen

- 1. Spülen Sie die Lösungsmittelbehälter und alle Teile der Lösungsmittelfilter gründlich aus. Verwenden Sie zum Spülen hochreines Lösungsmittel.
- 2. Setzen Sie den Lösungsmittelfilter zusammen; tragen Sie dabei geeignete saubere Handschuhe:
 - a) Setzen Sie den Schlauchadapter in den Filterhalter (Oberteil) ein.
 - b) Drehen Sie den Filterhalter (Oberteil) mit Adapter auf den Kopf und drücken Sie ihn gegen eine geeignete Oberfläche, so dass der Adapter fest im Filterhalter sitzt.
 - c) Legen Sie die Fritte waagrecht in den Filterhalter (Unterteil).
 - d) Schrauben Sie das Oberteil des Filters fest auf das Unterteil.



Abbildung 78: Lösungsmittelfilter

Nr.	Beschreibung
1	Schlauchadapter
2	Filterhalter, Oberteil
3	Filterfritte
4	Filterhalter, Unterteil

Gehen Sie wie folgt vor



Abbildung 79: Vorbereiten des Lösungsmittelschlauchs und des Behälterdeckels

Nr.	Beschreibung
1	Lösungsmittelschlauch
2	Schlauchführung
3	Behälterdeckel
4	Verschlusskappen
5	Lösungsmittelfilter mit Schlauchadapter

- 1. Führen Sie jeden Lösungsmittelschlauch durch eine Schlauchführung.
- 2. Führen Sie den Lösungsmittelschlauch durch eine Öffnung in einem Deckel für einen Lösungsmittelbehälter. Die Schlauchführung verhindert, dass der Lösungsmittelschlauch im Behälter verrutscht.
- 3. Verschließen Sie die Öffnungen in den Behälterdeckeln mit den Verschlusskappen.
- 4. Schieben Sie einen Lösungsmittelfilter auf jeden Lösungsmittelschlauch.
- 5. Füllen Sie die Lösungsmittelbehälter mit Lösungsmittel.
- Drehen Sie den Flaschendeckel handfest an. Drücken Sie die Schlauchführung in die Öffnung im Behälterdeckel, um den Schlauch im Deckel zu arretieren.
- 7. Stellen Sie die Lösungsmittelbehälter in das Solvent Rack.
- 8. Platzieren Sie die Lösungsmittelschläuche gerade in den Schlauchführungen.
- 9. Prüfen Sie die Lösungsmittelschläuche über den gesamten Flussweg hinweg. Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.

Sehen Sie dazu auch

Informationen zu Lösungsmitteln und Additiven (▶ Seite 144)

5.13 Füllen der Behälter für die Nadel-Waschflüssigkeit

Füllen Sie vor dem Betrieb des Systems die Behälter für die Nadel-Waschflüssigkeit.

TIPP Um die Bildung von kleinen Luftblasen in den Schläuchen zu vermeiden, platzieren Sie die Behälter im Solvent Rack oder oberhalb des Systemturms.

Erforderliche Werkzeuge

Kapillarschneider

Vorbereitungen

Spülen Sie die Behälter für die Waschflüssigkeit gründlich aus. Verwenden Sie zum Spülen hochreines Lösungsmittel.

Waschflüssigkeiten

Wählen Sie Waschflüssigkeiten, die für Ihre Anwendung geeignet sind und Probenreste zuverlässig von der Nadel entfernen. Es ist jeweils ein Behälter für starke und ein Behälter für schwache Waschflüssigkeit vorhanden. Typischerweise entspricht die starke Waschflüssigkeit dem Lösungsmittel B Ihrer Applikation oder ist ein stärkeres Lösungsmittel, während die schwache Waschflüssigkeit dem Anfangseluenten Ihres Gradienten ähnlich ist.

Art der Waschflüs sigkeit	Beschreibung (für Umkehrphasen- Anwendungen)	Schläuche
Stark (strong)	Waschflüssigkeit mit hohem organischen Anteil, zum Beispiel 0,1% Ameisensäure in Acetonitril	Waschschlauch für das Purgen und kurzzeitige Außerbetriebnahme mit Beschriftung S (rot): angeschlossen an Degaser-Eingang Strong in der Pumpe.
		Waschschlauch für das Waschen des Nadelinneren und -äußeren mit Beschriftung SWP (rot): angeschlossen an oberen Einlass der Nadel- Waschpumpe.
Schwach (weak)	Waschflüssigkeit mit niedrigem organischen Anteil, zum Beispiel 0,1% Ameisensäure in Wasser	Waschschlauch für das Waschen des Nadelinneren und das Purgen mit Beschriftung W (grün): angeschlossen an Degaser-Eingang Weak in der Pumpe.
		Waschschlauch für das Waschen des Nadeläußeren mit Beschriftung WWP (grün): angeschlossen an unteren Einlass der Nadel-Waschpumpe.

Gehen Sie für jeden Behälter wie folgt vor



Abbildung 80: Vorbereiten des Waschschlauchs und des Behälterdeckels

Nr.	Beschreibung
1	Waschschlauch
2	Schlauchführung
3	Behälterdeckel
4	Verschlusskappen
5	Lösungsmittelfilter mit Schlauchadapter

- 1. Führen Sie jeden Waschschlauch durch eine Schlauchführung.
- 2. Führen Sie jeden Waschschlauch durch eine Öffnung im Deckel eines der Behälter. Die Schlauchführung verhindert, dass der Lösungsmittelschlauch im Behälter verrutscht.
- 3. Verschließen Sie die Öffnungen in den Behälterdeckeln mit den Verschlusskappen.
- 4. Schieben Sie einen Filterhalter mit Adapter auf jeden Waschschlauch.
- 5. Befüllen Sie den Behälter mit Spülflüssigkeit.
- Drehen Sie den Flaschendeckel handfest an. Drücken Sie die Schlauchführung in die Öffnung im Behälterdeckel, um den Schlauch im Deckel zu arretieren.
- 7. Stellen Sie den Behälter in das Solvent Rack.
- 8. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.

Spülen des Nadel-Waschsystems

Die Nadel-Waschsysteme werden gespült, wenn Sie nach dem Einschalten des Systems das Skript **A02 - Autostart** ausführen. Siehe Vorbereiten des Systems für den ersten Betrieb (> Seite 138).

5.14 Füllen des Behälters für die Hinterspülflüssigkeit

Füllen Sie vor dem Betrieb des Systems den Behälter für die Hinterspülflüssigkeit.

TIPP Um die Bildung von kleinen Luftblasen in den Schläuchen zu vermeiden, platzieren Sie die Behälter im Solvent Rack oder oberhalb des Systemturms.

Erforderliche Werkzeuge

Kapillarschneider

Vorbereitungen

Spülen Sie den Hinterspülbehälter gründlich aus. Verwenden Sie zum Spülen hochreines Lösungsmittel.

Hinterspülflüssigkeit

Verwenden Sie 75% Isopropanol in Wasser und 0,1% Ameisensäure (in HPLC-Qualität; Aufbereitung nach Volumen, zum Beispiel 75 mL Isopropanol + 25 mL Wasser + 0,1 mL Ameisensäure).

Gehen Sie wie folgt vor





Nr.	Beschreibung
1	Spülschlauch
2	Schlauchführung
3	Behälterdeckel
4	Verschlusskappen

- 1. Führen Sie den Schlauch durch die Schlauchführung.
- 2. Führen Sie den Schlauch durch eine Öffnung im Deckel des Behälters. Die Schlauchführung verhindert, dass der Schlauch im Behälter verrutscht.
- 3. Verschließen Sie die Öffnungen im Behälterdeckel mit Verschlusskappen.
- 4. Befüllen Sie den Behälter mit Spülflüssigkeit.
- Drehen Sie den Flaschendeckel handfest an. Drücken Sie die Schlauchführung in die Öffnung im Behälterdeckel, um den Schlauch im Deckel zu arretieren.
- 6. Stellen Sie den Behälter in das Solvent Rack.
- 7. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.

5.15 Einschalten des Systems nach der Installation

Bevor Sie das System einschalten, installieren Sie den Säulenthermostat, insofern Ihre Systemkonfiguration einen Säulenthermostaten enthält.

Ein-/Ausschalter und Hauptnetzschalter

Die Abbildung zeigt die Ein-/Ausschaltelemente des Systems.



Abbildung 82: Ein-/Ausschalter und Hauptnetzschalter

Nr.	Beschreibung
1	System-Ein-/Ausschalter
2	Ein- und Ausschalter des System-Controllers
3	Hauptnetzschalter

Vorbereitungen

1. Stellen Sie sicher, dass die Transportsicherungsschraube des Autosamplers korrekt gelöst ist.

ACHTUNG—Schäden am System

Bevor Sie das Vanquish-System einschalten, vergewissern Sie sich, dass die Transportsicherungsschraube korrekt gelöst wurde, um Schäden am System zu vermeiden.

 Vergewissern Sie sich, dass die Chromatographie-Software auf dem Datensystemrechner installiert ist. Die erforderlichen USB-Treiber werden automatisch geladen und das Windows[™]-Betriebssystem kann die Systemmodule erkennen, wenn das System eingeschaltet wird.

Gehen Sie wie folgt vor

 Schalten Sie jedes Systemmodul (Pumpe, Autosampler, optionaler Säulenthermostat) über dessen Hauptnetzschalter ein. Schalten Sie den Autosampler ein, ehe Sie den Pumpenfluss einschalten und sich Druck aufbaut.

Die Hinterspülung wird über die Pumpe eingeschaltet und gesteuert. Jedes Mal, wenn Sie die Pumpe einschalten, führt sie zunächst einen Hinterspülzyklus durch.

- Schalten Sie den System-Controller über den Hauptnetzschalter ein. Um den System Controller einzuschalten, genügt es, den Hauptnetzschalter zu betätigen (Betätigung des Ein-/Ausschalters ist nicht nötig). Nachdem der System-Controller gebootet hat, leuchtet die Status-LED.
- 3. Schalten Sie das System über den System-Ein-/Ausschalter am Systemsockel an.

5.16 Integrieren in ein Geräte-LAN

Um eine Kommunikationsverbindung zwischen dem System und einem Geräte-LAN herzustellen, das aus dem Chromatographie-Datensystem und einem anderen Gerät besteht, zum Beispiel einem Massenspektrometer, müssen Sie das System in das Geräte-LAN integrieren. Das Vanquish User Interface bietet zwei Konnektivitätsoptionen:

- Über eine benutzerdefinierte IP-Adresse Wird zum Beispiel benötigt, wenn ein Gerät innerhalb eines Geräte-LANs eine benutzerdefinierte IP-Adresse besitzt.
- Über das Dynamic Host Configuration Protocol (DHCP)

Voraussetzungen

- Sie haben den System Controller mit dem Datensystemrechner verbunden.
- Sie haben auf dem Datensystemrechner, auf dem die Chromatographie-Software installiert ist, ein Geräte-LAN eingerichtet.
- *Nur bei benutzerdefinierten IP-Adressen*: Sie kennen den IP-Adressbereich und die Netzmaske des Geräte-LANs.

Gehen Sie wie folgt vor

- 1. Öffnen Sie das Einstellungsfenster auf dem Display.
- 2. Wählen Sie den Bereich Konnektivität.
- 3. Wählen Sie eine der Konnektivitätsoptionen aus.
- 4. Wenn Sie Benutzerdefiniert ausgewählt haben:
 - a) Geben Sie die Netzmaske des Geräte-LANs ein.
 - b) Geben Sie eine einzigartige IP-Adresse des System Controllers innerhalb Ihres Geräte-LANs ein.
- 5. Wählen Sie Anwenden aus.

Sehen Sie dazu auch

Anschließen eines LAN-Kabels (Seite 115)

5.17 Vorbereiten des Systems für den ersten Betrieb

ACHTUNG—Partikel und Luftblasen können die Systemfunktionalität beeinträchtigen

Einige Komponenten im System sind nach der Lagerung oder bei Auslieferung mit Isopropanol gefüllt.

- Um Luftblasen (und Isopropanol) vor der Erstinbetriebnahme des Systems auszuspülen, führen Sie das Skript aus: A02 - Autostart.
- Um Isopropanol auszuspülen, verwenden Sie Lösungsmittel, die mit Isopropanol mischbar sind. Ist dies nicht der Fall, stellen Sie das Lösungsmittel schrittweise um.

Wann

Vor dem ersten Gebrauch des Systems

Benötigte Mengen an Lösungsmitteln und Hinterspülflüssigkeit

Benötigte Flüssigkeit	Verbrauch während Aus- führung des Skripts: A02 - Autostart	Verbrauch während der Analyse	Anforderungen an die Flüssig- keiten
Lösungsmittel	Max. 10 mL je Lösungsmittel- behälter	Abhängig von der Flussrate	Wenn Sie das System oder eine Komponente zum ersten Mal in Betrieb nehmen: Verwenden Sie Lösungsmittel, die mit Isopropanol mischbar sind. Ist dies nicht der Fall, stellen Sie das Lösungsmittel schrittweise um. In allen anderen Fällen: Siehe Informationen zu Lösungsmitteln und Additiven (▶ Seite 144).
Hinterspül- flüssigkeit	Max. 10 mL	Circa 250 mL für 14 Tage	Siehe Füllen des Behälters für die Hinterspülflüssigkeit (▶ Seite 133).

Benötigte Mengen an Nadel-Waschflüssigkeit

Verbrauch wäh- rend Ausführung des Skripts: A02 - Autostart	Verbrauch wä lyse	hrend der Ana-	Anforderungen an die Flüssigkeiten		
Max. 10 mL	Verbrauch pro Nadel-Wasch- flüssigkeitsbehälter und pro Injektion ¹		Siehe Füllen der Behälter für die Nadel-Waschflüssigkeit (▶ Seite 130).		
	SWP	500 μL			
	WWP	400 µL			
	W	250 μL			
¹ : Die Werte gelten für: 25-uL Probenschleife. Einstellungen im Chromatographie-					

Die Werte geiten für: 25-μL Probenschleife, Einstellungen im Chromatographie Datensystem: direkte Injektion (Probenschleife in Bypass-Position), Standard-Wascheinstellungen.

Anmerkung: Der Verbrauch hängt von folgenden Faktoren ab: Probenschleifenvolumen, Säulengröße, Waschdauer, Workflow und Waschgeschwindigkeit. Wenn Sie andere Einstellungen verwenden, können die Werte wesentlich höher liegen.

Gehen Sie wie folgt vor

- Überprüfen Sie den Füllstand aller Behälter, vor allem den im Behälter für die Nadel-Waschflüssigkeit. Stellen Sie sicher, dass die Mengen ausreichen für:
 - Skript A02 Autostart (siehe Tabelle oben)
 - Erste Analyse (siehe Tabelle oben und addieren Sie den Verbrauch, der von der Flussrate Ihrer Anwendung abhängt).
- 2. Schließen Sie die Türen aller Module im Vanquish Neo-System, soweit noch nicht geschehen.
- Um das Lösungsmittel A/B einzustellen, führen Sie das Skript aus: A01 - Typen der Pumpenlösungsmittel einstellen.
- Führen Sie das Skript mit oder ohne Diagnosefunktion (Dichtigkeitstest) aus: A02 - Autostart. Das Skript spült den Systemflusspfad bis zur Säulen-Einlasskapillare, testet den Systemrückdruck und führt optional einen Dichtigkeitstest durch.
- 5. Um den Trennsäulentyp einzustellen, führen Sie das Skript aus: **A03 - Trennsäulentyp festlegen**.

- 6. Um die Trennsäuleneigenschaften einzustellen, führen Sie das Skript aus: **A04 Trennsäulenspezifikationen festlegen**.
- Um die Säule zu konditionieren, führen Sie das Skript aus: B06 Säule(n) konditionieren.

6 Betrieb

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

6.1 Sicherheitshinweise zum Betrieb

Beachten Sie beim Betrieb des Systems folgende Sicherheitshinweise:



Beachten Sie alle Warnhinweise und Vorsichtsmaßnahmen in Sicherheitsmaßnahmen (▶ Seite 26).



WARNUNG—Teile in Bewegung

Teile im Autosampler sind in Bewegung, wenn der Autosampler eine Probenanalyse vorbereitet und durchführt. Diese Teile können eine Klemmgefahr darstellen, die Verletzungen verursachen kann.

- Halten Sie die Autosamplertüren geschlossen, wenn der Autosampler eine Probenanalyse vorbereitet und durchführt.
- Während dieser Phasen leuchtet die LED-Leiste am Autosampler blau.



WARNUNG—Brennbare und Gefährliche Dämpfe von verschütteten Proben

Brennbare oder gefährliche Dämpfe von verschütteten Proben können sich im Autosampler ansammeln. Dies birgt Gefahren für Gesundheit und Sicherheit.

- Achten Sie darauf, dass Probenhalter und Probenplatten korrekt in den Segmenten stehen.
- Wenn Sie Probenhalter verwenden, setzen Sie die Probenfläschchen ein, bevor Sie die Probenhalter in den Drehteller stellen.
- Wenn Proben im Autosampler verschüttet wurden, schalten Sie den Autosampler aus. Wischen Sie die verschüttete Probe auf und lassen Sie die Autosamplertür offen. Warten Sie eine angemessene Zeit lang, bis die Stelle der verschütteten Probe getrocknet ist und sich sämtliche Dämpfe aufgelöst haben, bevor Sie den Autosampler wieder in Betrieb nehmen.



VORSICHT—Heiße Oberflächen

Oberflächen im Inneren des Systems 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:

- Um zu vermeiden, dass durch eventuelle Undichtigkeiten oder ein Trockenlaufen der Pumpe Schäden entstehen, stellen Sie immer die untere Druckgrenze für die Pumpe ein.
- Wenn eine Undichtigkeit im Gerät auftritt, schalten Sie den Pumpenfluss ab und beheben Sie umgehend die Ursache für die Undichtigkeit.
- Überzeugen Sie sich immer, dass der Autosampler eingeschaltet ist, ehe der Pumpenfluss an ist und sich Druck aufbaut. Ist der Autosampler ausgeschaltet, zum Beispiel aufgrund eines Stromausfalls, stoppen Sie den Pumpenfluss und warten Sie, bis der Druck auf null ist, bevor Sie den Autosampler und die anderen Module wieder einschalten.

6.2 Informationen zu Lösungsmitteln und Additiven

6.2.1 Allgemeine Kompatibilität

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

- Verwenden Sie nur Lösungsmittel und Additive, die mit allen Teilen im Flussweg kompatibel sind. Informationen zu den Materialien, die im Flussweg verwendet werden, finden Sie unter Leistungsspezifikationen (▶ Seite 390) für das jeweilige Modul.
- Das Material des Einsatzes f
 ür Flaschen besteht aus geschlossenzelligem expandiertem Polypropylen, das mit den g
 ängigsten HPLC-L
 ösungsmitteln kompatibel ist.
- Wasserfreies Methanol kann bei Titan-Oberflächen zu Spannungsbrüchen führen, insbesondere wenn dem Gemisch Ameisensäure oder TFA hinzugefügt wird. Thermo Fisher Scientific empfiehlt, 3% Wasser hinzuzufügen, um dem Problem vorzubeugen.

Kompatibilität der Kolbendichtungen

- Bei Auslieferung der Pumpe sind Dichtungen aus UHMW-PE installiert. Bei Verwendung von Tetrahydrofuran, Ketonen oder Ammoniumhydroxid als Lösungsmittel können die Dichtungen beschädigt werden.
- 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.

6.2.2 Erlaubte pH-Bereiche

Erlaubter pH-Bereich für das System:

Erlaubter pH-Bereich	Bemerkung
2-10	<i>pH-Werte von 2</i> : Die Anwendungsdauer sollte möglichst kurz sein. Spülen Sie das System nach entsprechenden Anwendungen gründlich.
6.2.3 Erlaubte Konzentrationen

Konzentration	Wert	
Pufferkonzentration	k.A.	
Chloridkonzentration:	Weniger oder gleich 1 mol/L Hohe Chlorid-Konzentration: Die Anwendungsdauer sollte möglichst kurz sein. Spülen Sie das System nach entsprechenden Anwendungen gründlich.	

Erlaubte Konzentrationen für das System:

6.2.4 Verwenden von Lösungsmitteln und Additiven

Partikel, die in das Chromatographie-System gelangen, können Kapillaren und Ventile blockieren, den Verschleiß erhöhen und die Säule oder das System beschädigen. Speziell in wässrigen Lösungsmitteln können sich Algen und andere Mikroorganismen vermehren, sich im System ablagern und die Lösungsmittelfilter verstopfen. Verstopfte Kapillaren und Filter können zu erhöhtem oder instabilem Systemdruck führen.

Beachten Sie im Hinblick auf die optimale Leistungsfähigkeit des Systems die folgenden Hinweise.

Lösemittelqualität und Filtration

- Verwenden Sie je nach Erfordernis Ihrer Anwendung hochreine Lösungsmittel (gefiltert) und Additive, zum Beispiel in UHPLC-Qualität oder LC/MS-Qualität. Gefilterte hochreine Lösungsmittel sind von den Herstellern in der Regel entsprechend gekennzeichnet.
- Verwenden Sie hochwertiges Wasser, zum Beispiel in UHPLC-Qualität oder LC/MS-Qualität (0,2 μm gefiltert).
- Bei der Verwendung von Wasser aus Wasseraufbereitungsanlagen kann es zu polymeren Verunreinigungen kommen, wenn die Aufbereitungsanlage nicht ordnungsgemäß gewartet ist.
- Wenn Sie Salzlösungen oder Puffer vorbereiten und die Vorbereitung abgeschlossen ist, verwenden Sie Membranfiltration (0,2 μm), um Schmutzpartikel zu entfernen und mikrobielles Wachstum zu reduzieren.
- Verwenden Sie die von Thermo Fisher Scientific empfohlenen Lösungsmittelfilter. Prüfen Sie die Filterfritten regelmäßig auf Durchlässigkeit und tauschen Sie die Fritten gegebenenfalls aus.

- Verwenden Sie frisches, gut entgastes Lösungsmittel. Tauschen Sie die Lösungsmittel mindestens wöchentlich aus. Tauschen Sie das Lösungsmittel im Behälter stets vollständig aus. Füllen Sie keine Lösungsmittelreste auf, um Schwankungen in der Lösungsmittelzusammensetzung oder die Bildung ungewollter Bestandteile zu vermeiden.
- Recyceln Sie im Hinblick auf die längere Lebensdauer der Dichtungen keine Lösungsmittel und fördern Sie nicht im Kreislauf.
- Beachten Sie die spezifischen Eigenschaften der Lösungsmittel, wie Viskosität, Mischbarkeit, Siedepunkt, UV-Absorption, sowie die Löslichkeit von Salzen in jedem Verhältnis im Verlauf des Gradienten.

Lösungsmittelbehälter

- Bevor Sie einen Lösungsmittelbehälter befüllen, spülen Sie den Behälter gründlich mit einem hochreinen Lösungsmittel aus.
- Zur Verringerung von Algenwachstum verwenden Sie braune Glasbehälter. Verwenden Sie geeignete Additive für wässrige Lösungsmittel, zum Beispiel Ameisensäure.
- Vermeiden Sie die Verwendung von Reinigungsmittel f
 ür die Reinigung von Glasbeh
 ältern. Sp
 ülen Sie Glasbeh
 älter f
 ür LC/MS-Anwendungen (einschlie
 ßlich Messzylinder) vor der Verwendung mit L
 ösungsmitteln in LC/MS-Qualit
 ät; zudem sollten diese separat beschriftet und gelagert werden.
- Verwenden Sie kein Methanol aus Aluminiumbehältern.

Spülen nach dem Betrieb

- Spülen Sie Puffer und peroxid-bildende Lösungsmittel aus.
- Lassen Sie keine Puffer, Salzlösungen oder aggressive Lösungsmittel ohne Fluss über einen längeren Zeitraum im System stehen.
- Bevor Sie das Laufmittel von einem Puffer oder einer Salzlösung auf ein organisches Lösungsmittel umstellen, spülen Sie das System gründlich mit Wasser in UHPLC-Qualität oder LC/MS-Qualität durch.
- Bevor Sie auf einen anderen Lösungsmitteltyp umstellen, vergewissern Sie sich, dass das neue Lösungsmittel mit dem vorherigen Lösungsmittel mischbar ist. Sind die Laufmittel nicht mischbar, kann es zu Ausflockungen kommen.

 Wechseln Sie zwischen zwei nicht miteinander mischbaren Lösungsmitteln mit Hilfe eines Löslichkeitsvermittlers, zum Beispiel Isopropanol. Ersetzen Sie das erste Lösungsmittel schrittweise durch den Löslichkeitsvermittler, und ersetzen Sie dann schrittweise den Löslichkeitsvermittler durch das neue Lösungsmittel.

6.3 Allgemeine Hinweise zum Systembetrieb

Beachten Sie die folgenden Informationen und Hinweise zum Betrieb des Systems.

So verbessern Sie die Leistung des Systems

- Waschen Sie die Nadel von innen und außen, um Verschleppung zu reduzieren. Das Waschen der Nadel stellt sicher, dass die Nadel ordnungsgemäß von jeglichen Probenresten gereinigt wird.
- Halten Sie während des Betriebs die Tür des Säulenthermostaten geschlossen.
- Vermeiden Sie das Trockenschalten der Ventile bei abgeschaltetem Pumpenfluss. Dadurch kann das Ventil beschädigt werden.
- Betreiben Sie die Pumpe nicht mit einer trockenen Hinterspülung.
- Beachten Sie die allgemeinen Richtlinien und Empfehlungen zur Verwendung von Lösungsmitteln und Additiven im System. Siehe Informationen zu Lösungsmitteln und Additiven (▶ Seite 144).

Hinweise zur Probenbearbeitung

Beachten Sie beim Betrieb des Autosamplers Folgendes:

TIPP—Autosamplertüren während der Probenbearbeitung geschlossen halten

- Wenn die Türen während der Bewegung der Nadel und des Drehtellers geöffnet werden, hält der Autosampler die laufende Bewegung des Nadelarms an. Die Nadel sticht nicht in das Probengefäß ein und zieht keine Probe auf. Es ertönt ein akustisches Signal, das anzeigt, dass die Türen offen sind.
- Jedes Mal, wenn die T
 üren geschlossen werden, f
 ührt der Autosampler einen 'Inventory Scan' der Probenhalter und -gef
 äße im Probenraum durch.

Sehen Sie dazu auch

Optimieren der Leistung (> Seite 167)

6.4 Einschalten des Systems

- Schalten Sie den System-Controller über den oberen Ein-/ Ausschalter am Systemsockel ein. Nachdem der System-Controller gebootet hat, leuchtet die Status-LED.
- 2. Schalten Sie das System über den unteren System-Ein-/Ausschalter am Systemsockel an.

6.5 Bedienelemente

Das System wird hauptsächlich über einen Rechner gesteuert, auf dem die Chromatographie-Software installiert ist, sowie über das Vanquish User Interface.

Zusätzlich stehen folgende Bedienelemente am System zur Verfügung:

- Tastaturen: Über die Tasten auf den Tastaturen können Sie einige Funktionen direkt am System ausführen.
- Statusanzeigen:
 Über die Statusanzeigen können Sie schnell den Betriebszustand des Gerätes erkennen. Im System gibt es folgende Statusanzeigen:
 - LED-Statusleiste vorne unten an jedem Gerät.
 - **STATUS**-LEDs auf den Tastaturen.
 - Status-LED am System-Controller

6.5.1 Tastaturen

Die folgenden Elemente stehen zur Verfügung:

• STATUS-LED

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.

Tasten

Über die Tasten können Sie einige Funktionen direkt am Gerät ausführen. Beim Drücken einer Taste bestätigt ein akustisches Signal, dass die Aktion ausgeführt wird.

STATUS-LED

Die **STATUS**-LED auf der Tastatur 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 die Meldung auf dem Benachrichtigungsfenster. Abhilfemaßnahmen finden Sie unter Fehlersuche (> Seite 343).	

Informationen zu den LEDs neben einer Taste auf der Tastatur finden Sie in der Tabelle zu den Tasten.

Tasten auf der Tastatur

Die Tasten haben folgende Funktion:

Autosampler und Pumpe

Taste	Funktion		
DOCK	Durch Drücken dieser Taste wird der gewählte Pumpenkopf in die geeignete Position für Wartungsarbeiten gebracht, zum Beispiel für den Tausch von Pumpenkopf oder Kolben.		
	Wenn sich der Pumpenkopf in der Position für Wartungsarbeiten befindet, beginnt die LED neben der Taste FLOW grün zu blinken. Durch erneutes Drücken der Taste wird der Pumpenkopf in die Betriebsposition zurückgefahren. Die LED blinkt nicht mehr grün.		
FLOW	 Durch Drücken der Taste wird der Fluss gestartet oder gestoppt, wobei die gewählte Flussrate und die Lösungsmittelzusammensetzung berücksichtigt werden. Die LEDs L und R neben der Taste beziehen sich auf den linken bzw. rechten Pumpenkopf: LED aus = Pumpenfluss ist ausgeschaltet oder Null LED grün = Pumpenkopf fördert LED blinkt grün = Pumpenkopf befindet sich in der Position für Wartungsarbeiten 		
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 innerhalb von 10 Minuten. Andernfalls ertönt das Signal erneut. Wenn das Gerät ein anderes Problem erkennt, ertönt sofort wieder ein akustisches Signal.		
PURGE	Reserviert für zukünftige Verwendung. Informationen zum Purgen finden Sie unter Purgen (Übersicht) (▷ Seite 65).		
ROTATE	Durch Drücken der Taste können Sie den Drehteller gegen den Uhrzeigersinn zur nächsten Beladungsposition drehen.		
SELECT	Autosampler • Reserviert für zukünftige Verwendung. <i>Pumpe</i> Durch Drücken der Taste wird der Pumpenkopf ausgewählt, für den Sie eine DOCK -Funktion über die Tastatur ausführen können:		
	 LED L grün = linker Pumpenkopf ausgewählt LED R grün = rechter Pumpenkopf ausgewählt 		
SERVICE	 Die Taste hat folgende Funktionen: Durch einmaliges Drücken der Taste fährt die Probennadel vom Nadelsitz hoch und das Injektionsventil schaltet in die Position Bypass. Dies ist zum Beispiel erforderlich, um die Nadeleinheit durch Festziehen der Transportsicherungsschraube zu sichern. Wenn Sie die Taste ein zweites Mal drücken, wird die Nadel initialisiert und die Nadel bewegt sich über den Waschport. Wenn Sie die Taste ein drittes Mal drücken, fährt die Nadel zurück in 		
	den Nadelsitz und die Nadel wird initialisiert. Hinweis: Die Transportsicherungsschraube muss gelöst sein.		

Taste	Funktion
VALVE	Durch Drücken der Taste können Sie das Injektionsventil in die Position Bypass schalten. Die LED B leuchtet grün, wenn sich das Injektionsventil in der Position Bypass befindet.
WASH	Das Drücken der Taste startet einen Waschzyklus des Nadeläußeren, in dem die Nadel von außen im Waschport gewaschen wird. Nach dem Nadel-Waschzyklus fährt die Nadel zurück zum Nadelsitz.

Säulenthermostat

Taste	Funktion		
MUTE ALARM	Ein akustisches Signal ertönt, wenn das Gerät ein Problem erkannt hat. Drücken Sie diese Taste, um das akustische Signal für den aktuellen Alarm auszuschalten. Beseitigen Sie die Ursache innerhalb von 10 Minuten. Andernfalls ertönt das Signal erneut. Wenn das Gerät ein anderes Problem erkennt, ertönt sofort wieder ein akustisches Signal.		
TEMP CTRL	 Durch Drücken der Taste wird die Temperaturregelung ein- oder ausgeschaltet: LED neben der Taste leuchtet grün = Temperaturregelung ist eingeschaltet LED neben der Taste aus = Temperaturregelung ist ausgeschaltet 		
VALVE LEFT	Durch Drücken der Taste schalten Sie das linke Säulenschaltventil (falls ein Ventil installiert ist). Durch Drücken und Halten der Taste starten Sie die Funktion "Catch Valve", wenn ein VICI-Ventil in der linken Ventilposition installiert ist. Bei 2-Positions 6-Port-Ventilen geben die LEDs neben den Tasten Auskunft über die Ventilposition:		
	 Obere LED grün = Ventil ist in Position 1 Untere LED grün = Ventil ist in Position 2 LEDs aus = Ventil schaltet 		
VALVE RIGHT	 Durch Drücken der Taste schalten Sie das rechte Säulenschaltventil (falls ein Ventil installiert ist). Durch Drücken und Halten der Taste starten Sie die Funktion "Catch Valve", wenn ein VICI-Ventil in der rechten Ventilposition installiert ist. Bei 2-Positions 6-Port-Ventilen geben die LEDs neben den Tasten Auskunft über die Ventilposition: Obere LED grün = Ventil ist in Position 1 Untere LED grün = Ventil ist in Position 2 LEDs aus = Ventil schaltet 		

6.5.2 LED-Statusleisten

Bei geschlossenen Gerätetüren geben die LED-Statusleisten an der Vorderseite des Moduls Auskunft. Ist das System im Chromatographie-Datensystem verbunden, zeigt die LED-Leiste gegebenenfalls weniger Informationen an.

Autosampler

LED-Leiste	Beschreibung		
Aus (dunkel)	Das Modul ist ausgeschaltet.		
Gedimmt	Die Modultüren sind geöffnet.		
Gelb, langsam blinkend	Das Modul ist eingeschaltet oder äquilibriert, aber das System ist nicht im Chromatographie-Datensystem verbunden.		
Gelb	Das System ist im Chromatographie-Datensystem verbunden, aber das Modul ist nicht äquilibriert.		
Grün, blinkend	Das Modul äquilibriert. Wenn Sie die Probenraum-Thermostatisierung nutzen, zeigt diese Farbe an, dass die Temperatur für die Thermostatisierung noch nicht erreicht ist.		
Grün	Das System ist im Chromatographie-Datensystem verbunden und das Modul ist äquilibriert, aber es läuft keine Datenaufnahme. Wenn Sie die Probenraum-Thermostatisierung nutzen, zeigt diese Farbe an, dass die Temperatur für die Thermostatisierung erreicht ist.		
Blau, Lauflicht	Der Autosampler führt Vorbereitungen für die Injektion durch. Teile des Autosamplers sind in Bewegung.		
Blau	Das System ist im Chromatographie-Datensystem verbunden und eine Injektion oder Sequenz läuft, einschließlich Datenaufnahme. Oder: Ein Diagnosetest läuft.		
Rot	Ein Problem oder Fehler ist aufgetreten. Prüfen Sie die Meldung auf dem Benachrichtigungsfenster. Abhilfemaßnahmen finden Sie unter Fehlersuche (> Seite 343).		

Pumpe

LED-Leiste	Beschreibung		
Aus (dunkel)	Das Modul ist ausgeschaltet.		
Gedimmt	Die Modultüren sind geöffnet.		
Gelb, langsam blinkend	Das Modul ist eingeschaltet oder äquilibriert, aber das System ist nicht im Chromatographie-Datensystem verbunden.		
Gelb	Das System ist im Chromatographie-Datensystem verbunden, aber das Modul ist nicht äquilibriert. Der Pumpenfluss ist aus.		
Grün, blinkend	Es läuft ein Spülzyklus.		
Grün	Das System ist im Chromatographie-Datensystem verbunden und das Modul ist äquilibriert, aber es läuft keine Datenaufnahme. Der Pumpenfluss ist an.		
Blau	Das System ist im Chromatographie-Datensystem verbunden und eine Injektion oder Sequenz läuft, einschließlich Datenaufnahme. Oder: Ein Diagnosetest läuft.		
Rot	Ein Problem oder Fehler ist aufgetreten. Prüfen Sie die Meldung auf dem Benachrichtigungsfenster. Abhilfemaßnahmen finden Sie unter Fehlersuche () Seite 343).		

Säulenthermostat

LED-Leiste	Beschreibung	
Aus (dunkel)	Das Modul ist ausgeschaltet.	
Gedimmt	Die Modultüren sind geöffnet.	
Gelb, langsam blinkend	Das Modul ist eingeschaltet oder äquilibriert, aber das System ist nicht im Chromatographie-Datensystem verbunden.	
Gelb	Das System ist im Chromatographie-Datensystem verbunden, aber das Modul ist nicht äquilibriert.	
Grün, blinkend	Das Modul äquilibriert.	
Grün	Das System ist im Chromatographie-Datensystem verbunden und das Modul ist äquilibriert, aber es läuft keine Datenaufnahme.	
Blau	Das System ist im Chromatographie-Datensystem verbunden und eine Injektion oder Sequenz läuft, einschließlich Datenaufnahme. Oder: Ein Diagnosetest läuft.	
Rot	Ein Problem oder Fehler ist aufgetreten. Prüfen Sie die Meldung auf dem Benachrichtigungsfenster. Abhilfemaßnahmen finden Sie unter Fehlersuche (> Seite 343).	

6.6 Vanquish User Interface

6.6.1 System Controller-Informationen

Die System Controller-Informationen werden in der oberen linken Ecke angezeigt.



Abbildung 83: System Controller-Informationen

Nr.	Beschreibung	
1	Uhrzeit ¹	
2	Datum ¹	
3	Firmware-Version des System Controllers	
4	Seriennummer des System Controllers	
¹ : Die Darstellung hängt von der Spracheinstellung ab.		

6.6.2 Systemstatus-Symbole

Die folgenden Symbole zeigen den Systemstatus an.

Symbol	Systemstatus
	ОК
	Fehler Ein oder mehrere Systemmodule haben einen Fehler ausgegeben und die LED-Leiste eines oder mehrerer Systemmodule leuchtet rot.
	Informationen zur Behebung des Fehlerzustands finden Sie unter System-Fehlersuche (> Seite 369).

6.6.3 Navigationsleistensymbole und andere Symbole

Die folgenden Symbole erscheinen in der Navigationsleiste.

Symbol	Beschreibung
	Startfenster Erlaubt den Zugriff auf die Funktionsparameter aller Systemmodule, die über die System-Interlink-Schnittstelle an den System Controller angeschlossen sind. Weitere Informationen finden Sie unter Startfenster (> Seite 156).
×	Skriptfenster Beinhaltet automatisierte Skripte. Weitere Informationen finden Sie unter Skriptfenster () Seite 157).
Ļ	Benachrichtigungsfenster Speichert alle Warnungen und Fehler der Systemmodule und des System Controllers seit dem letzten Booten des System Controllers Weitere Informationen finden Sie unter Meldungen (> Seite 346).
\$ 0	Einstellungsfenster Weitere Informationen finden Sie unter Einstellungsfenster (▶ Seite 157).
()	Informationsfenster Zeigt verschiedene Informationen an, zum Beispiel, alle angeschlossenen Systemmodule und ihre Seriennummern
M	Online-Plot-Fenster (verfügbar im Startfenster) Öffnet ein Dialogfeld, das die verbundenen Module in der oberen Hälfte und die dazugehörigen physikalischen Werte in der unteren Hälfte anzeigt. Das Diagramm wird kontinuierlich aktualisiert und zeigt das Signal der letzten 10 Minuten an.
∇	Filter (nur verfügbar im Skriptfenster und im Benachrichtigungsfenster) Ermöglicht zum Beispiel das Filtern nach einem Systemmodul.

6.6.4 Betriebszustand des Systems

Wenn das System im Chromatographie-Datensystem verbunden ist, wird der Betriebszustand (zum Beispiel Leerlauf oder Waschen) des Systems in der oberen rechten Ecke des Bildschirms angezeigt.

6.6.5 Startfenster

Im Startfenster wird jedes Systemmodul durch einen Bereich mit einem spezifischen Symbol dargestellt, welches folgende Informationen enthält:

• Funktionsparameter, zum Beispiel den Pumpendruck

 Jeder Bereich hat am unteren Rand einen bunten Balken, der die LED-Leiste darstellt. Eine Beschreibung der LED-Leistenfarben finden Sie in LED-Statusleisten (> Seite 153).



Abbildung 84: Systemmodul-Bereich (hier: Pumpe)

6.6.6 Skriptfenster

Das Skriptfenster stellt automatisierte Skripte für verschiedene Produktlebensphasen zur Verfügung, zum Beispiel für die Inbetriebnahme und die Fehlersuche.

Wenn Sie ein Skript anwählen, öffnet sich ein Dialogfenster, in dem Sie detaillierte Informationen und Anweisungen erhalten.

6.6.7 Einstellungsfenster

In diesem Fenster können Sie Folgendes tun:

- Vanquish User Interface konfigurieren:
 - Sprache
 - Einheiten
 - Zeitzone
 - Konnektivitätsmodus
- Ansehen:
 - Systemname, Firmware-Version und Betriebsstundenzahlen
 - Software-Lizenzen

6.7 Anpassen des Displays

6.7.1 Mögliche Positionen des Displays

Das Display kann auf verschiedene Weise positioniert werden:

- Vordere Position
 - Nach hinten gekippt



6.7.2 Verstellen des Displays von der vorderen Position in die seitliche Position

1. Drehen sie das Display um 180 Grad nach links.



Abbildung 85: Drehen des Displays nach links um 180 Grad

2. Klappen Sie das Display nach unten.



Abbildung 86: Klappen des Displays nach unten

6.8 Vorbereiten des Systems für den Betrieb

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

- 1. Spülen Sie das System. Siehe Spülen des Systems (> Seite 160).
- Äquilibrieren Sie das System. Siehe Äquilibrieren des Systems (> Seite 162).
- Beladen Sie den Drehteller. Siehe Beladen des Drehtellers (▶ Seite 163).
- Beachten Sie die Empfehlungen zu Kappen und Verschlüssen von Probengefäßen. Siehe Empfehlungen zu Kappen und Verschlüssen von Probengefäßen (▶ Seite 166).

6.8.1 Spülen des Systems

ACHTUNG—Partikel und Luftblasen können die Systemfunktionalität beeinträchtigen

Einige Komponenten im System sind nach der Lagerung mit Isopropanol gefüllt.

- Um Luftblasen (und Isopropanol) vor der erneuten Inbetriebnahme des Systems oder nach der Installation von zusätzlichen Komponenten in den Flusspfad auszuspülen, führen Sie das Skript aus: A02 - Autostart.
- Um Isopropanol auszuspülen, verwenden Sie Lösungsmittel, die mit Isopropanol mischbar sind. Ist dies nicht der Fall, stellen Sie das Lösungsmittel schrittweise um.

Wann

- Nach dem Einschalten des Systems
- Nach der Installation von Komponenten in den Flusspfad

Benötigte Flüssigkeit	Verbrauch während Aus- führung des Skripts: A02 - Autostart	Verbrauch während der Analyse	Anforderungen an die Flüssig- keiten
Lösungsmittel	Max. 10 mL je Lösungsmittel- behälter	Abhängig von der Flussrate	Wenn Sie eine Komponente zum ersten Mal in Betrieb nehmen: Verwenden Sie Lösungsmittel, die mit Isopropanol mischbar sind. Ist dies nicht der Fall, stellen Sie das Lösungsmittel schrittweise um. In allen anderen Fällen: Siehe Informationen zu Lösungsmitteln und Additiven (▶ Seite 144).
Hinterspül- flüssigkeit	Max. 10 mL	Circa 250 mL für 14 Tage	Siehe Füllen des Behälters für die Hinterspülflüssigkeit (▶ Seite 133).

Benötigte Mengen an Lösungsmitteln und Hinterspülflüssigkeit

Benötigte Mengen an Nadel-Waschflüssigkeit

Verbrauch wäh- rend Ausführung des Skripts: A02 - Autostart	Verbrauch während der Ana- lyse		Anforderungen an die Flüssigkeiten
Max. 10 mL	Verbrauch pro Nadel-Wasch- flüssigkeitsbehälter und pro Injektion ¹ :		Siehe Füllen der Behälter für die Nadel-Waschflüssigkeit (▶ Seite 130).
	SWP	500 μL	
	WWP	400 μL	
	W	250 μL	

¹: Die Werte gelten für: 25-µL Probenschleife, Einstellungen im Chromatographie-Datensystem: direkte Injektion (Probenschleife in Bypass-Position), Standard-Wascheinstellungen.

Anmerkung: Der Verbrauch hängt von folgenden Faktoren ab:

Probenschleifenvolumen, Säulengröße, Waschdauer, Workflow und Waschgeschwindigkeit. Wenn Sie andere Einstellungen verwenden, können die Werte wesentlich höher liegen.

Gehen Sie wie folgt vor

- Überprüfen Sie den Füllstand aller Behälter, vor allem den im Behälter für die Nadel-Waschflüssigkeit. Stellen Sie sicher, dass die Mengen ausreichen für:
 - Skript A02 Autostart (siehe Tabelle oben)
 - Erste Analyse (siehe Tabelle oben und addieren Sie den Verbrauch, der von der Flussrate Ihrer Anwendung abhängt).
- 2. Schließen Sie die Türen aller Module im Vanquish Neo-System, soweit noch nicht geschehen.
- Wenn Sie den Lösungsmitteltyp ändern möchten: Um den Lösungsmitteltyp A/B einzustellen, führen Sie das Skript aus: A01 -Typen der Pumpenlösungsmittel einstellen.
- 4. Dieser Schritt ist abhängig von der Situation:
 - Wenn Sie die zuvor installierte Säule verwenden: Pr
 üfen Sie die Säulen-Konfiguration und passen Sie diese gegebenenfalls mit den Skripten an: A03 - Trennsäulentyp festlegen und A04 -Trennsäulenspezifikationen festlegen.
 - Wenn Sie eine neue, unbenutzte Säule installiert haben: Konfigurieren Sie die neue Säule mit den Skripten: A03 -Trennsäulentyp festlegen und A04 -Trennsäulenspezifikationen festlegen.
- Führen Sie das Skript mit oder ohne Diagnosefunktion (Dichtigkeitstest) aus: A02 - Autostart. Das Skript spült den Systemflusspfad bis zur Säulen-Einlasskapillare, testet den Systemrückdruck und führt optional einen Dichtigkeitstest durch.

6.8.2 Äquilibrieren des Systems

Wann

- Nach Einschalten des Systems und nach dem Vorbereiten des Systems f
 ür die Inbetriebnahme
- Nach einer kurzzeitigen Außerbetriebnahme
- Nach Änderung der Flussrate
- Nach Änderung der Temperatur-Einstellungen

Gehen Sie wie folgt vor

Die System-Äquilibrierung sollte Folgendes umfassen:

1. Heizen (oder Kühlen) aller temperaturgeregelten Geräte für eine bestimmte Zeit:

Äquilibrierungsschritte	Dauer
Heizen (oder Kühlen) aller temperaturgeregelten Geräte im System auf die Anfangstemperatur. Temperaturgeregelte Geräte können zum Beispiel sein:	5 Stunden, abhängig von den Temperaturunterschieden, die für die Anwendung benötigt werden
 Pumpen-Flowmeter 	
 Thermostatisierter Probenraum im Autosampler 	
 Säulenthermostat und Säule 	

- 2. Gehen Sie während des Heizens (oder Kühlens) abhängig von der Situation wie folgt vor:
 - Wenn Sie die zuvor installierte Säule verwenden: Um die Säule zu äquilibrieren, führen Sie das Skript aus: B02 - Säule(n) reinigen und äquilibrieren.
 - Wenn Sie eine neue, unbenutzte Säule installiert haben: Um neue Säulen zu konditionieren, führen Sie das Skript aus: B06 – Säule(n) konditionieren.
- Beobachten Sie die Drift des Pumpendrucks und überprüfen Sie, dass die Drift 5-24 Stunden lang stabil in einem vernünftigen Bereich für die Anwendung, abhängig von der Retentionszeitgenauigkeit, die für die Anwendung benötigt wird.

6.8.3 Beladen des Drehtellers

Der Drehteller kann mit bis zu vier Probenhaltern und/oder Mikrotiterplatten bestückt werden.

Erforderliche Teile

Probenfläschchen, Probenhalter bzw. Mikrotiterplatten

Beachten Sie die Empfehlungen unter Empfehlungen zu Kappen und Verschlüssen von Probengefäßen (▶ Seite 166).

Vorbereitungen

- 1. Positionieren Sie die Fläschchen mit Ihren Proben im Probenhalter.
- 2. Überprüfen Sie die eingestellte Temperatur für den Probenraum und passen Sie sie gegebenenfalls an. Siehe Thermostatisierung des Autosampler-Probenraums (▶ Seite 173).
- Prüfen Sie die LED-Statusleiste am Autosampler. Öffnen Sie die Gerätetüren nicht, wenn die LED-Leiste blau leuchtet. Teile des Autosamplers sind in Bewegung.

Gehen Sie wie folgt vor

- 1. Drehen Sie das gewünschte Segment des Drehtellers nach vorn. Sie können den Drehteller folgendermaßen drehen:
 - Drücken Sie die Taste ROTATE auf der Tastatur. Dadurch wird der Drehteller gegen den Uhrzeigersinn zur nächsten Beladungsposition gedreht.
 - Drehen Sie den Drehteller manuell in die gewünschte Richtung.
 - Drehen Sie den Drehteller vom Vanquish Neo-ePanel im Chromatographie-Datensystem oder vom Autosampler-Bereich auf dem Startfenster im Vanquish User Interface aus.
- Stellen Sie einen Probenhalter oder eine Mikrotiterplatte in das ausgewählte Segment. Der Probenhalter bzw. die Mikrotiterplatte muss im Ausrichtungsrahmen des Segments sitzen und auf den Ausrichtungspunkten aufliegen.

ACHTUNG—Korrekte Ausrichtung von Probenhaltern

Werden Probenhalter und Mikrotiterplatten falsch ausgerichtet, kann dies zu Beschädigungen am Autosampler oder zu falschen Ergebnissen führen.

Stellen Sie Probenhalter und Mikrotiterplatten immer in richtiger Ausrichtung in den Drehteller, sodass sich Position A1 vorne links am Probenhalter bzw. der Platte befindet.



Abbildung 87: Korrekte Ausrichtung im Drehteller (hier mit einem Probenhalter)

- 3. Bestücken Sie die Zusatz-Fläschchenhalter, falls erforderlich.
- 4. Drehen Sie den Drehteller zur nächsten Beladungsposition und beladen Sie den Drehteller wie gewünscht. Beladen Sie auf diese Weise alle weiteren Segmente nach Erfordernis.
- Schließen Sie die Autosamplertüren. Der Autosampler beginnt einen 'Inventory Scan' der Probenhalter und -gefäße im Probenraum.

Automatische Probenhalter-Format-Erkennung

Wenn Sie Probenhalter oder Mikrotiterplatten mit Barcode zur Vanquish-Probenhalter-Format-Erkennung verwenden, werden das Format und die Ausrichtung des Probenhalters bei der Bestandsprüfung automatisch erkannt.

Probenhalter-Format manuell setzen

Wenn Sie Probenhalter oder Mikrotiterplatten ohne speziellen Vanquish-Barcode verwenden, müssen Sie das Probenhalter-Format für jedes Segment auf dem Vanquish Neo-ePanel im Chromatographie-Datensystem einstellen.

Sehen Sie dazu auch

- B Drehteller (► Seite 50)
- Probenhalterformat-Erkennung (> Seite 51)

6.8.4 Empfehlungen zu Kappen und Verschlüssen von Probengefäßen

Verwenden Sie ausschließlich die hier beschriebenen Kappen für Probenfläschchen und Abdeckungen für Mikrotiterplatten.

Kappen und Septen für Probenfläschchen

Verwenden Sie nur talkumfreie ungefüllte Silikon/PTFE-basierte Septen. Talkum und Partikel aus Dichtungen können Probleme mit Rückdruck erzeugen oder das System verstopfen.

Beispiel für Kappen, die die Anforderungen erfüllen: Thermo Scientific Schraubkappen und Septen mit rotem PTFE oder weißem Silikon

Verschlussmatten für Mikrotiterplatten

Verwenden Sie ungefüllte Silikon-Verschlussmatten.

Beispiel für Matten, die die Anforderungen erfüllen: Thermo Scientific WebSeal[™] 96-Well unbeschichtete Mikrotiterplatten aus Kunststoff mit WebSeal Silikonmatten

Dichtungsbänder für Mikrotiterplatten

Verwenden Sie Dichtungsbänder bzw. Wiederversiegelungsbänder ohne Klebstoff im Kavitätenbereich. Klebstoff im Kavitätenbereich kann zu einer Kontamination der Injektionsnadel führen.

Beispiel für Bänder, die die Anforderungen erfüllen: Thermo Scientific WebSeal 96-Well unbeschichtete Mikrotiterplatten aus Kunststoff mit WebSeal Dichtungsband mit 3 nicht-klebenden Schichten im Kavitätenbereich

Verwenden Sie *keine* Versiegelungen, bei denen die Gefahr besteht, dass Dichtungsmaterial durch die Probennadel ausgestochen wird. Zum Beispiel können Aluminiumteilchen aus Aluminium-Versiegelungsfolien ausgestochen werden und sich im Nadelsitz ablagern.

6.9 Optimieren der Leistung

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

6.9.1 Abhängigkeit von verfügbarem Pumpendruck und Flussrate

Der maximale Druck, den die Pumpe liefern kann, hängt wie folgt ab:

- Eingestellte Flussrate
- Viskosität des Lösungsmittels, das in einem der beiden Kanäle verwendet wird (Kanal A oder Kanal B)

Für die vorkalibrierten Lösungsmittel zeigt der Graph, wie der maximal verfügbare Pumpendruck von der eingestellten Flussrate und dem Kanal abhängt.



Abbildung 88: Abhängigkeit von verfügbarem Pumpendruck und Flussrate für Kanal A und Kanal B

Beachten Sie für andere Lösungsmittel Folgendes:

- Verwenden Sie f
 ür Gradienten-Flussraten bis 10 μL/min, das wässrige (schwache) L
 ösungsmittel in Kanal A und das organische (starke) L
 ösungsmittel in Kanal B.
- Wenn die Gesamtflussrate 10 μL/min übersteigt und der erwartete Systemrückdruck 140 MPa (1400 bar) übersteigt, verwenden Sie das Lösungsmittel mit der höchsten Viskosität in Kanal A.

Im Betrieb überprüft die Pumpe kontinuierlich, ob sie die eingestellte Flussrate bei dem sich daraus ergebenden Druck liefern kann. Wenn der Systemrückdruck so hoch wird, dass die Pumpe die eingestellte Flussrate nicht liefern kann, passiert Folgendes:

- Die Injektion oder Sequenz wird abgebrochen.
- Es erscheint eine Meldung, die zum Beispiel darauf hinweist, dass die Pumpe den geforderten Fluss bei diesem Druck nicht liefern konnte (Code 4531) oder dass der Druck den absoluten Grenzwert überschritten hat (Code 4503).

Wenn Sie die Flussrate reduzieren, nimmt der Systemrückdruck ab. Gleichzeitig nimmt der von der Pumpe verfügbare Druck zu. Wenn die Pumpe die eingestellte Flussrate immer noch nicht liefern kann, reduzieren Sie den Systemrückdruck weiter. Zum Beispiel, indem Sie die Flussrate weiter reduzieren, die Säulentemperatur erhöhen oder den Innendurchmesser der Flusskapillaren erhöhen.

6.9.2 Hinzufügen von kundenspezifischen Lösungsmitteln zur Kalibriertabelle

Die Pumpe kann exakte Flussraten nur liefern, wenn das Lösungsmittel, das Sie verwenden möchten, kalibriert wurde. Kalibrierte Lösungsmittel können im Skript **A01 - Typen der Pumpenlösungsmittel einstellen** oder in der Kalibriertabelle der Lösungsmittel ausgewählt werden. Für die gängigsten Lösungsmitteltypen und Gemische wurde die Kalibrierung im Werk durchgeführt:

- Wasser, 100%
- Acetonitril, 100%
- Gemisch aus Acetonitril und Wasser, 80:20 (v:v)
- Gemisch aus Methanol und Wasser, 90:10 (v:v)

Wenn Sie andere Lösungsmittel verwenden möchten, kalibrieren Sie diese einmal, um sie der Konfiguration hinzuzufügen. Die Kalibrierung wird für einen Lösungsmitteltyp empfohlen, der > 5% (volumenbezogen) von einem der oben genannten Typen abweicht.

Beispiele

Beispiel 1

Lösungsmittel		Vorgehen	
A	98:2 (v:v), Wasser:Acetonitril plus 0,1% Ameisensäure	Verwenden Sie die Kalibrierdaten von 100% Wasser.	
В	80:20 (v:v), Acetonitril:Wasser plus 0,1% Ameisensäure	Verwenden Sie die Kalibrierdaten von 80:20% Acetonitril.	

Beispiel 2

Lösungsmittel		Vorgehen	
A	95:5 (v:v), Wasser:Dimethylsulfoxid plus 0,1% Ameisensäure	Verwenden Sie die Kalibrierdaten von 100% Wasser.	
В	75:20:5 (v:v:v), Acetonitril:Wasser:Dimethylsulfoxid plus 0,1% Ameisensäure	Verwenden Sie die Kalibrierdaten von 80:20% Acetonitril.	

Beispiel 3

Lösungsmittel		Vorgehen	
A	100% Wasser plus 0,1% Ameisensäure	Verwenden Sie die Kalibrierdaten von 100% Wasser.	
В	90:10 (v:v), Acetonitril:Wasser plus 0,1% Ameisensäure	Kalibrierung ist erforderlich. Siehe weiter unten.	

Allgemeine Hinweise

- Die Lösungsmittelkalibrierung erfolgt pro Kanal.
- Wenn Sie ein Lösungsmittel in beiden Kanälen verwenden möchten, müssen Sie die Kalibrierung zweimal durchführen, um das Lösungsmittel für jeden Kanal separat zu kalibrieren.
- Die eigentliche Kalibrierung ist automatisiert und läuft unbeaufsichtigt. Bevor Sie das Skript für die eigentliche Kalibrierung starten, müssen Sie die Pumpe für die Kalibrierung einrichten und konfigurieren und Benutzereingaben vornehmen. Zum Abschluss müssen Sie die ursprüngliche Kapillarkonfiguration der Pumpe wiederherstellen und alle Kapillaren und das System wieder entsprechend der Anwendung verbinden.
- Kalibrierskripte führen Sie durch alle Schritte des Vorgangs.
- Der eigentliche Kalibriervorgang dauert circa 100 Minuten. Planen Sie außerdem im Vorfeld 30 Minuten für die Vorbereitung des Kalibrieraufbaus ein und 30 Minuten danach für die Wiederherstellung der Betriebskonfiguration.

Erforderliche Teile und Utensilien

- Lösungsmittel, geeignet zum Purgen, um schädliche Substanzen herauszuspülen
- Lösungsmittel, das Sie kalibrieren möchten
- Viper-Verschlussfittinge (3 Fittinge)

Gehen Sie wie folgt vor

- 1. Setzen Sie auf dem **Skriptfenster** den Filter auf **Lösungsmittelkalibrierung**.
- 2. Führen Sie die Skripte für die Lösungsmittelkalibrierung nacheinander aus und folgen Sie den Anweisungen.
- Nachdem Sie die Betriebskonfiguration wieder hergestellt haben, testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: E04 -Detaillierter Dichtigkeitstest (Pumpe).

Ergebnisse der Kalibrierung

- Kalibrierung erfolgreich: Das Lösungsmittel wird in der Kalibriertabelle hinterlegt und kann ausgewählt werden.
- Während der Kalibrierung ist ein Fehler aufgetreten: Schließen Sie über die Kalibrierskripte entweder die Kalibrierung ab oder stellen Sie die Betriebskonfiguration wieder her.

6.9.3 Ändern des Lösungsmitteltyps

- 1. Wenn Sie den Lösungsmitteltyp ändern möchten:
 - a) Um den Lösungsmitteltyp A/B einzustellen, führen Sie das Skript aus: **A01 Typen der Pumpenlösungsmittel einstellen**.
 - b) Um das System zu spülen, führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit den folgenden Parametern: Lösungsmittel A/B und Nur aktualisieren = Aus).

6.9.4 Ändern der Hinterspülflüssigkeit

Wenn Sie die Hinterspülflüssigkeit ändern möchten:

Führen Sie das Skript aus: **B01 - Flüssigkeiten/Lösungsmittel ändern** (mit den folgenden Parametern: **Hinterspülflüssigkeit hinten** und **Nur aktualisieren = Aus**).

6.9.5 Ändern der Nadel-Waschflüssigkeit

Wenn Sie die Nadel-Waschflüssigkeit ändern möchten:

Führen Sie das Skript aus: **B01 - Flüssigkeiten/Lösungsmittel ändern** mit den entsprechenden Nadel-Waschparametern und **Nur aktualisieren = Aus**.

6.9.6 Durchführen des Nullpunktabgleichs für die Flusssensoren der Pumpe

Das Skript zur Durchführung des Nullpunktabgleichs der Flusssensoren misst das Sensorsignal bei einem Fluss von Null und speichert das Ergebnis, damit das Flusssignal entsprechend korrigiert werden kann. Damit wird die Nullpunktverschiebung, die seit der Kalibrierung im Werk aufgetreten ist, ausgeglichen.

Abgleich nicht erforderlich

Ein Nullpunktabgleich der Flusssensoren muss nicht durchgeführt werden, wenn die Retentionszeiten stabil sind oder im erwarteten Bereich liegen. Führen Sie den Nullpunktabgleich der Flusssensoren in diesen Situationen nicht durch:

- Während eine wissenschaftliche Studie oder Messreihe läuft.
- Während einer Messreihe, in der vergleichbare Ergebnisse erwartet werden.
- Sehr häufig, zum Beispiel, täglich.

Abgleich kann erforderlich sein

Im Hinblick auf genaue und vergleichbare Messergebnisse kann es in diesen Situationen erforderlich sein, einen Nullpunktabgleich der Flusssensoren durchzuführen:

- Wenn die Retentionszeiten deutlich vom erwarteten Bereich abweichen (verschoben sind). Dies kann dazu führen, dass Peaks fehlen.
- Wenn Retentionszeiten im Vergleich zu früheren Messungen stark variieren.

Der Nullpunktabgleich der Flusssensoren muss schätzungsweise einmal alle 1 bis 3 Monate durchgeführt werden, besonders bei Analysen < 2 µL/min.

Gehen Sie wie folgt vor

Beachten Sie Folgendes, ehe Sie beginnen:

- Das System muss mindestens 5 Stunden eingeschaltet sein. Nur dann ist der Abgleich möglich.
- *Empfehlung*: Führen Sie den Abgleich mit den Lösungsmitteln Ihrer Applikation durch.
- Führen Sie für den Abgleich dieses Skript aus: C01 -Nullpunktabgleich des Pumpenflusssensors durchführen.
- 2. Planen Sie nach dem Abschluss des Abgleichs etwas Äquilibrierzeit für eine optimale Retentionszeitstabilität ein, zum Beispiel eine Blindprobe (blank run).

6.9.7 Waschen der Probennadel

Der Nadelwaschzyklus wäscht die Nadel, um Probenreste von der Nadel zu entfernen. Dies kann Teil der Probensequenz sein oder manuell durchgeführt werden. Dieser Abschnitt beschreibt, wie Sie einen Nadel-Waschzyklus manuell auslösen, und damit die Systemleistung und die Workflow-Abfolge optimieren.

Wann

- Austauschen der Nadel-Waschflüssigkeit
- Tauschen der Waschschläuche
- Regelmäßige und vorbeugende Wartungsarbeiten für das Nadel-Waschsystem

Vorbereitungen

Beachten Sie die Hinweise zum Waschen der Nadel unter Füllen der Behälter für die Nadel-Waschflüssigkeit (▶ Seite 130).

Gehen Sie wie folgt vor

So starten Sie das Waschen des Nadeläußeren:

Drücken Sie die Taste **WASH** auf der Tastatur.

Dies startet einen Waschzyklus des Nadeläußeren, in dem die Nadel von außen im Waschport gewaschen wird.

So starten Sie das Waschen des Nadelinneren und -äußeren manuell:

Führen Sie auf dem Vanquish User Interface folgendes Skript aus: **B03 -** System reinigen

Das System wäscht daraufhin das Äußere sowie das Innere der Nadel.

Nachdem der Nadel-Waschzyklus abgeschlossen wurde, fährt die Nadel zurück zum Nadelsitz.

Sehen Sie dazu auch

- Nadel-Waschsystem (Übersicht) (▶ Seite 58)
- Wartung des Nadel-Waschsystems (> Seite 270)

6.9.8 Thermostatisierung des Autosampler-Probenraums

Im Autosampler kann der Probenraum thermostatisiert werden.

Gehen Sie wie folgt vor

- 1. Aktivieren Sie die Temperaturregelung im Chromatographie-Datensystem oder im Autosampler-Bereich auf dem Startfenster im Vanquish User Interface.
- 2. Legen Sie die Solltemperatur fest. Prüfen Sie, dass die Thermostatisierungstemperatur für Ihre Proben geeignet ist.

TIPP Der Autosampler speichert beim Ausschalten die Solltemperatur. Überprüfen Sie die Temperatur beim Einschalten des Autosamplers und passen Sie sie gegebenenfalls an.

6.9.9 Thermostatisierung des Säulenraums

Der Säulenraum kann thermostatisiert werden.

Gehen Sie wie folgt vor

- 1. Schalten Sie die Temperaturregelung im Vanquish User Interface oder durch Drücken der Taste **TEMP CTRL** ein.
- 2. Wählen Sie die Solltemperatur im Chromatographie-Datensystem oder im Säulenthermostat-Bereich auf dem Startfenster im Vanquish User Interface. Prüfen Sie, dass die Thermostatisierungstemperatur für folgende Teile geeignet ist:
 - Ventile
 - Säulen

Hinweise zum Einstellen der Temperatur

Der Säulenthermostat hat kein aktives Kühlsystem. Beachten Sie Folgendes:

- Sie können die Temperatur im Säulenthermostaten nicht unter die Temperatur senken, die sich aus der Umgebungstemperatur und der Eigenerwärmung des Gerätes ergibt.
- Der Temperatursollwert muss immer mindestens 5 °C höher als die Umgebungstemperatur eingestellt werden.

Der Zustand Ready		
	Das Property Ready im Chromatographie-Datensystem gibt an, ob der Säulenthermostat betriebsbereit ist (Ready). Solange der Säulenthermostat nicht betriebsbereit ist (NotReady), können Sie nicht mit der Analyse beginnen. Die Einstellungen, die Sie für Ready Temp Delta und Equilibration Time vornehmen, beeinflussen die Zeit, die der Säulenthermostat benötigt, um in den Zustand Ready zu gelangen.	
Ready Temp Delta		
	Ready Temp Delta gibt an, um wieviel Grad die aktuelle Temperatur über bzw. unter der vorgegebenen Solltemperatur liegen darf. Weicht die aktuelle Temperatur um mehr als den hier eingegebenen Wert von der Solltemperatur ab, geht der Säulenthermostat in den Zustand NotReady über und ist nicht betriebsbereit.	
	Beispiel: Temperatursollwert: 45 °C, Ready Temp Delta: 1 °C	
	Der Säulenthermostat ist betriebsbereit (Ready), wenn die aktuelle Temperatur für die Equilibration Time zwischen 44 °C und 46 °C liegt. Liegt die aktuelle Temperatur außerhalb dieses Bereichs, geht der Säulenthermostat in den Zustand NotReady über.	
Äquilibrierdauer		
	Gibt an, wie lange die gewünschte Temperatur, unter Berücksichtigung der Einstellung unter Ready Temp Delta , stabil bleiben muss, ehe der Säulenthermostat betriebsbereit ist und mit der Analyse begonnen werden kann.	
	Beispiel: Temperatursollwert: 45 °C, Ready Temp Delta : 1 °C, Equilibration Time : 0,5 min	
	Der Säulenthermostat ist betriebsbereit (Ready), wenn die aktuelle Temperatur zwischen 44 °C und 46 °C liegt und 0,5 Minuten gehalten wird.	
	Beachten Sie folgende Hinweise:	
	• Je kleiner der gewählte Wert für Ready Temp Delta und je größer der Wert für Equilibration Time ist, desto länger dauert es, bis der Säulenthermostat betriebsbereit ist und Sie mit der Datenaufnahme bzw. Analyse beginnen können.	
	 Äußere Umstände, wie Zugluft oder Klimaanlagen, können dazu führen, dass der Säulenthermostat gar nicht oder nur nach sehr langer Äquilibrierzeit in den Zustand Ready übergeht. 	
	• Beachten Sie, je besser das System äquilibriert ist, desto besser sind die chromatographischen Ergebnisse.	

6.10 Außerbetriebnehmen des Systems

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

6.10.1 Kurze Betriebsunterbrechung des Systems

Beachten Sie folgende Hinweise für das System, wenn Sie den Betrieb für kurze Zeit, zum Beispiel über Nacht, unterbrechen (kurzzeitige Außerbetriebnahme):

Vorbereitungen

Setzen Sie die Maximaltemperatur auf 40 °C, insofern Ihre Systemkonfiguration einen Säulenthermostaten enthält.

Gehen Sie wie folgt vor

1. Führen Sie das Skript aus und folgen Sie den Anweisungen: **B05 -**Außerbetriebnahme (kurzzeitig).

Inbetriebnehmen des Systems nach der kurzzeitigen Außerbetriebnahme

- 1. Gehen Sie wie folgt vor, bevor Sie den Betrieb wieder aufnehmen:
 - a) Lassen Sie Pumpenfluss und Temperatur auf die Betriebsbedingungen zurück äquilibrieren.
 - b) Spülen Sie den Autosampler mit schwacher Nadelwaschflüssigkeit. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern mit den Parametern Waschen des Nadelinneren, schwache Flüssigkeit und Nur aktualisieren = An.
- 2. Führen Sie das Skript aus: **B02 Säule(n) reinigen und äquilibrieren**.

6.10.2 Lange Betriebsunterbrechung des Systems

Folgen Sie bei längeren Betriebsunterbrechungen, bei denen das System im Labor verbleibt, den folgenden Anweisungen.

1. Führen Sie das Skript aus: **B05 - Außerbetriebnahme (langfristig)** mit dem folgenden Parameter: **Zu entfernendes Modul = Keines**.

TIPP Die Lösungsmittelschläuche müssen in der Regel nicht aus den Schlauchführungen entfernt werden (auch nicht, wenn der Moduleinschub getauscht wird). Wenn Sie jedoch die Lösungsmittelschläuche aus den Schlauchführungen in der Pumpe und allen Modulen oberhalb der Pumpe im Systemturm entfernen, achten Sie darauf, nicht an den anderen Schläuchen in den Führungen zu ziehen.

- 2. Leeren Sie die Lösungsmittelbehälter.
- 3. Schalten Sie das System über den System-Ein-/Ausschalter (unterer Schalter am Systemsockel) aus.
- 4. Schalten Sie den System-Controller über seinen Ein-/Ausschalter (oberer Schalter am Systemsockel) aus.

TIPP Wenn der Autosampler ausgeschaltet wird, öffnet sich die linke Tür des Autosamplers automatisch, um eine ordnungsgemäße Belüftung des Probenraums zu ermöglichen. Die Tür lässt sich nicht schließen, solange der Autosampler ausgeschaltet ist.

Inbetriebnehmen des Systems nach der langfristigen Außerbetriebnahme

- 1. Füllen Sie die Lösungsmittelbehälter. Siehe Füllen der Lösungsmittelbehälter (▶ Seite 128).
- Schalten Sie das System ein. Siehe Einschalten des Systems nach der Installation (▶ Seite 135).
- 3. Bereiten Sie das System für den Betrieb vor. Siehe Vorbereiten des Systems für den Betrieb (▶ Seite 160).

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).

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 325).

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 26).



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—Scharfe Spitze der Injektionsnadel

Die Injektionsnadel besitzt eine scharfe Spitze, die Verletzungen der Haut verursachen kann.

Berühren Sie niemals die Nadelspitze, um Personenschäden zu vermeiden.



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 des Systems 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. Siehe Entleeren der Lösungsmittelschläuche (> Seite 234).
- 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.
ACHTUNG

Selbst wenn Sie den Vanquish-Systemturm nur geringfügig auf dem Arbeitstisch verschieben, kann der Systemsockel beschädigt und dessen Funktionalität beeinträchtigt werden.

Bevor Sie den Systemturm verschieben, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln.

TIPP Wenn der Autosampler ausgeschaltet wird, öffnet sich die linke Tür des Autosamplers automatisch, um eine ordnungsgemäße Belüftung des Probenraums zu ermöglichen. Die Tür lässt sich nicht schließen, solange der Autosampler ausgeschaltet ist.

7.3 Allgemeine Regeln für Wartung und Service

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

Allgemeine Regeln

- 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.

Öffnen von Flussverbindungen

- 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.
- Nachdem Sie Wartungs- oder Servicearbeiten an den Flussverbindungen im System vorgenommen haben, gehen Sie wie folgt vor: Testen Sie zunächst den Systemrückdruck (Skript: D01 -Systemrückdruck testen). Testen Sie dann das System auf Undichtigkeiten (Skript: D02 - Systemdichtigkeit testen (was getestet werden soll = System)).

ACHTUNG

Flussverbindungen und Kapillaranschlüsse sind äußerst empfindlich gegen Verunreinigungen. Staub und Verschmutzungen können diese Anschlüsse verunreinigen.

Verschließen Sie Kapillaren stets mit Schutzkappen und offene Flussverbindungen mit Verschlussstopfen, um sie vor Verunreinigungen zu schützen.

Reparatur

• Wenden Sie sich an den Thermo Fisher Scientific-Kundendienst vor Ort, wenn Sie das System für eine Reparatur an Thermo Fisher Scientific zurückschicken müssen.

Sehen Sie dazu auch

Ersatzteile und Verbrauchsmaterialien (> Seite 406)

7.4 Wartung und Wartungsintervalle

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

7.4.1 Wartungszeitplan

Führen Sie regelmäßig die Wartungsarbeiten in den folgenden Tabellen durch. Die in den Tabellen 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 System verwendet werden.

7.4.1.1 Wöchentliche Wartungsarbeiten

Produktkom- ponente	Was ist zu tun
System/Alle Module	 Flussverbindungen: Prüfen Sie die Flussverbindungen auf Anzeichen von:
	 Blockierung: Führen Sie das Skript aus und folgen Sie den Anweisungen: D01 - Systemrückdruck testen.
	 Salzablagerungen
	 Luftblasen Entgasen Sie die Lösungsmittel mit Ultraschall und purgen Sie die Pumpe und den Autosampler, wenn nötig.
	 Lösungsmittel: Wenn Sie eines der folgenden Lösungsmittel verwenden, gehen Sie folgendermaßen vor:
	 Reines Wasser oder ein wässriger Puffer: Um den Lösungs- mittelbehälter auszutauschen, führen Sie dieses Skript aus: B01 Flüssigkeiten/Lösungsmittel ändern.
	 Puffer oder Salzlösungen: Spülen Sie das System nach Ab- schluss der Arbeiten gründlich mit einem geeigneten Lösungs- mittel, das keine Puffer oder Salze enthält. Siehe Purgen (Übersicht) (> Seite 65).
	 Korrosive Lösungsmittel oder Methanol mit <10% Wasser: Spülen Sie das System mit Wasser oder mit anderen geeigneten, nichtkorrosiven Lösungsmitteln. Siehe Purgen (Übersicht) (> Seite 65).

Produktkom- ponente	Was ist zu tun
Waschsysteme	 Überprüfen Sie den Füllstand. Überprüfen Sie, ob sich Schmutzpartikel, Staub oder Algen in den Waschbehältern befinden. Wenn sich Schmutzpartikel, Staub oder Algen in den Waschbehältern befinden, ersetzen Sie die Waschbehälter durch saubere Waschbehälter. Führen Sie das Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern (Nur aktualisieren = An). Stellen Sie sicher, dass die Waschbehälter frei von Reinigungsmittel- oder Chemikalienrückständen sind. Wir empfehlen, die Waschbehälter mit einer kleinen Menge der geplanten Waschflüssigkeit zu spülen.
Autosampler	 Prüfen Sie die Probenfläschchen und Mikrotiterplatten im Autosampler auf Risse oder Schäden. Wischen Sie verschüttete Proben im Autosampler auf, falls erforderlich.

7.4.1.2 Regelmäßige Wartungsarbeiten

Produktkom- ponente	Was ist zu tun
System/Alle Module	 Tauschen Sie die Lösungsmittelbehälter alle 1-2 Wochen aus. Um Lösungsmittelbehälter auszutauschen, führen Sie dieses Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern. Spülen Sie die Gefäße vor dem Auffüllen gründlich aus.
	 Prüfen Sie das System auf Undichtigkeit: D02 - Systemdichtigkeit testen (was getestet werden soll = System).
	 Erneuern Sie die Filterfritten der Lösungsmittelschläuche in regelmäßigen Abständen. Dies ist speziell bei der Verwendung von wässrigen Lösungsmitteln wichtig. In wässrigen Lösungs- mitteln können sich mit der Zeit Algen und andere Mikro- organismen vermehren und die Filterfritten verstopfen.
	 Pr üfen Sie die Flussverbindungen auf Anzeichen von Besch ädigungen, wie Knicke, Risse, Schnitte oder Blockierung.
	 Pr üfen Sie die Systemdrainage auf Blockierungen. Siehe Testen des Drainagesystems (
	Leeren Sie die Abfallbehälter.
	 Reinigen Sie das System. Siehe Reinigen oder Dekontaminieren eines Systemmoduls oder des Systems () Seite 190).
	 Pr üfen Sie die Warnaufkleber am System und an den System- modulen. Sollte ein Warnaufkleber fehlen oder nicht mehr deutlich lesbar sein, wenden Sie sich f ür Ersatzaufkleber an Thermo Fisher Scientific.

Produktkom- ponente	Was ist zu tun
Waschsysteme	 Nadel-Waschsysteme: Ersetzen Sie die Behälter für die Nadel-Waschflüssigkeit regel- mäßig, ungefähr alle 1-2 Wochen. Um die Behälter für die Nadel- Waschflüssigkeit zu tauschen, führen Sie dieses Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern.
	• Hinterspülung:
	 Tauschen Sie den Behälter für die Hinterspülflüssigkeit aus. Spülen Sie den Behälter gründlich, bevor Sie ihn befüllen. Verwenden Sie zum Spülen hochreines Lösungsmittel.
	 Prüfen Sie die Hinterspülschläuche. Wenn ein Schlauch blockiert (gequetscht oder verbogen) ist, tauschen Sie den Schlauch aus.
	 Überprüfen Sie den Peristaltikschlauch in der Hinterspülpumpe. Wenn er verrutscht oder verhärtet ist, korrigieren Sie die Position oder ersetzen Sie den Schlauch. Siehe Tauschen des Peristaltikschlauchs in der Hinterspülpumpe (* Seite 289). Tauschen Sie die Schläuche alle sechs Monate als präventive Wartungsmaßnahme aus.
	 Testen Sie die Hintersp ülung auf Dichtigkeit. Siehe Testen der Hintersp ülung auf Undichtigkeit (Seite 279).
Pumpe	 Reinigen Sie die Pumpenkopf-Kugelventile einmal pro Monat. Siehe Reinigen der Kugelventile (Pumpenkopf) (
	 Um eine Kontamination zu vermeiden, spülen Sie die Pumpe mit frischen Lösungsmitteln. Siehe Spülen der Pumpe zur Vermeidung von Kontamination (▶ Seite 187).
Autosampler	 Tauschen Sie den Drainagepumpen-Schlauch alle sechs Monate aus. Siehe Tauschen des Drainagepumpen-Schlauchs () Seite 267).
	 Nur für den Workflow vom Typ Trap und Elution (beheizt): Bestimmen Sie den Rückdruck des Trap-Säulenflusspfads. Führen Sie das Skript aus: D01 - Systemrückdruck testen mit dem Parameter: Nur Trap-Säulen-Widerstand testen. Notieren Sie diesen Wert und vergleichen Sie ihn mit nachfolgenden Rückdrucktests, um die Alterung der Trap-Säule zu beurteilen.
Display	Ziehen Sie die Schrauben nach, die das Display am System be- festigen.

Sehen Sie dazu auch

Verwenden von Lösungsmitteln und Additiven (> Seite 145)

7.4.1.3 Jährliche Wartungsarbeiten

Produktkom- ponente	Was ist zu tun
System/Alle Module	Lassen Sie einmal pro Jahr vorbeugende Wartungsmaßnahmen von einem Thermo Fisher Scientific-Servicetechniker durchführen.
Hinterspülung	 Tauschen Sie die Hinterspüldichtungen. Siehe Tauschen der Hinterspüldichtung (> Seite 215). Tauschen Sie die Hinterspülschläuche aus. Siehe Wartung der Hinterspülung (> Seite 279).
Pumpe	Prüfen Sie die Kolben optisch auf Ablagerung von Partikeln oder Anzeichen einer Beschädigung, zum Beispiel Kratzer oder Risse. Siehe Reinigen des Kolbens (▶ Seite 202).

7.4.2 Spülen der Pumpe zur Vermeidung von Kontamination

Erforderliche Teile

Wie zutreffend:

- Frisches Lösungsmittel
- Isopropanol
- Ethanol

Beachten Sie die folgenden Hinweise

Die folgenden Hinweise beziehen sich auf die konventionelle Verwendung des Vanquish-Systems wie in dieser Anleitung empfohlen, zum Beispiel unter Verwendung von hochwertigen Lösungsmitteln (mindestens UHPLC-Qualität oder LC/MS-Qualität), die häufig ausgetauscht werden:

- Im Normalfall ist es ausreichend, die Pumpe regelmäßig mit frischem Lösungsmittel zu spülen.
- Wenn Sie Wasser, Acetonitril oder Methanol f
 ür Ihre Chromatographie verwenden, sollten Sie die Pumpe ein Mal pro Woche sp
 ülen.
- Mobile Phasen, die Salze, Puffersubstanzen und/oder andere Additive enthalten, sind grundsätzlich anfällig für organische Verschmutzung und/oder biologischen Abbau. Wenn Sie nichtflüchtige Pufferlösungen verwenden, beachten Sie Folgendes:
 - Mobile Phasen müssen vor ihrer Verwendung gefiltert werden.
 - Für einen einwandfreien Betrieb muss das System wöchentlich gespült werden, zum Beispiel mit > 60% Isopropanol oder Ethanol.

• Passen Sie das Intervall an die verwendeten Lösungsmittel an.

TIPP Ziehen Sie bei hartnäckiger Verschmutzung, zum Beispiel bei reproduzierbar auftretenden Störpeaks im Chromatogramm ohne Injektion einer Probe, eine schärfere Reinigungsprozedur in Betracht. Siehe Spülen bei hartnäckiger Verschmutzung (> Seite 189).

Sehen Sie dazu auch

Verwenden von Lösungsmitteln und Additiven (> Seite 145)

7.4.3 Spülen bei hartnäckiger Verschmutzung

Wann

TIPP Bei konventioneller Verwendung des Vanquish Neo-Systems, wie in dieser Anleitung empfohlen, ist es im Normalfall ausreichend, die Pumpe regelmäßig zu spülen. Siehe Spülen der Pumpe zur Vermeidung von Kontamination () Seite 187).

Ziehen Sie zum Beispiel bei reproduzierbar auftretenden Störpeaks im Chromatogramm ohne Injektion einer Probe, eine schärfere Reinigungsprozedur in Betracht. Folgen Sie den Anweisungen.

Erforderliche Teile und zusätzliche Utensilien

- Geeignete Gegendruckkapillare, die circa 20 90 MPa bei einem Fluss von 50 μL/min mit Wasser erzeugt (zum Beispiel, nanoViper-Kapillare, Fused Silica, Innendurchmesser: 20 μm, Länge: 350 mm)
- 6 molare (6 M) Salpetersäure (HPLC-Qualität)
- Abfallbehälter (zum Auffangen der Salpetersäure)
- Frisches Wasser (HPLC-Qualität)
- Frisches Acetonitril (HPLC-Qualität)
- Neue Lösungsmittelbehälter mit frischem Lösungsmittel
- Neue Fritten für die Lösungsmittelfilter

Gehen Sie wie folgt vor

- 1. Installieren Sie die Gegendruckkapillare am Pumpenausgang. Führen Sie das offene Ende der Kapillare in den Abfallbehälter.
- 2. Tauschen Sie die Fritten in den Lösungsmittelfiltern gegen neue Fritten.
- 3. Tauschen Sie den installierten Abfallbehälter gegen den neuen Abfallbehälter aus, um chemische Reaktionen der Salpetersäure mit dem Lösungsmittel im Abfallbehälter zu vermeiden.
- 4. Spülen Sie die Pumpe mit 6-molarer Salpetersäure für 2 Stunden mit 50% B bei einer Flussrate von 50 μ L/min.
- 5. Spülen Sie die Pumpe mit frischem Wasser in HPLC-Qualität, bis der pH-Wert neutral ist.
- 6. Entfernen Sie den Abfallbehälter mit der Salpetersäure und installieren Sie wieder den zuvor installierten Abfallbehälter.

- 7. Spülen Sie die Pumpe mit frischem Acetonitril für 2 Stunden mit 50% B bei einer Flussrate von 50 μ L/min.
- 8. Schließen Sie die neuen Lösungsmittelbehälter an, die frisches Lösungsmittel enthalten.
- 9. Entfernen Sie die Gegendruckkapillare und schließen Sie das System wieder entsprechend Ihrer Anwendung an.
- 10. Purgen Sie die Pumpe.
- 11. Äquilibrieren Sie das System.

Sehen Sie dazu auch

Purgen (Übersicht) (Seite 65)

7.4.4 Reinigen oder Dekontaminieren eines Systemmoduls oder des Systems

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 eines Systemmoduls oder des Systems. Verwenden Sie ein geeignetes Reinigungs- oder Desinfektionsmittel, um sicherzustellen, dass die Systemmodule oder das System nach der Behandlung sicher gehandhabt werden können.

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 Gerätes gelangen. Flüssigkeiten können einen Kurzschluss auslösen, wenn sie in Kontakt mit elektronischen Komponenten kommen.

Vorbereitungen

- 1. Nur, wenn Sie den Autosampler oder das System reinigen oder dekontaminieren wollen:
 - a) Fahren Sie die Nadeleinheit in die Service-Position.
 - b) Ziehen Sie die Transportsicherungsschraube fest (siehe Sichern der Nadeleinheit (▶ Seite 248)).
 - c) Nehmen Sie alle Probengefäße und Probenhalter aus dem Probenraum.

- 2. Fahren Sie nach Erfordernis fort:
 - Wenn Sie ein einzelnes Systemmodul reinigen oder dekontaminieren wollen: Schalten Sie das Systemmodul aus und trennen Sie das Netzkabel von der Stromversorgung.
 - Wenn Sie das System reinigen oder dekontaminieren wollen: Schalten Sie alle Systemmodule und das System aus und trennen Sie die 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.

Schritte nach der Reinigung/Dekontamination

- 1. Nur, wenn Sie den Autosampler oder das System gereinigt oder dekontaminiert haben:
 - a) Lösen Sie die Transportsicherungsschraube.
 - b) Stellen Sie die Probengefäße und Probenhalter in den Probenraum.
- 2. Schließen Sie das/die Netzkabel an die Stromversorgung an.
- Schalten Sie das/die Systemmodul(e) ein.
 Wenn der Autosampler eingeschaltet ist, fährt die Nadeleinheit automatisch in die Position Inject.
- 4. Bereiten Sie das System für den Betrieb vor (siehe Vorbereiten des Systems für den Betrieb (▶ Seite 160)).

7.5 Wartung des Pumpenkopfs

Dieser Abschnitt beschreibt, wie der Pumpenkopf oder Pumpenkopfteile gewartet oder getauscht werden. Folgen Sie den Schritten für die Wartungsprozedur, die Sie ausführen möchten.

- Zum Tausch eines Pumpenkopfs siehe Tauschen des Pumpenkopfs (> Seite 195).
- Zum Tausch eines Kolbens siehe Tauschen der Kolben (> Seite 200).
- Zum Reinigen eines Kolbens siehe Reinigen des Kolbens (▶ Seite 202).
- Zum Tausch eines Stützrings siehe Tauschen des Stützrings (> Seite 203).
- Zum Tausch einer Kolbendichtung siehe Tauschen der Kolbendichtung (▶ Seite 208).
- Zum Tausch einer Hinterspüldichtung siehe Tauschen der Hinterspüldichtung (▶ Seite 215).
- Zum Tausch einer Pumpenkopfdichtung siehe Tauschen der Pumpenkopfdichtungen (▶ Seite 219).
- Zum Reinigen oder zum Tausch eines Kugelventils siehe Warten der Pumpenkopf-Kugelventile (▶ Seite 224).

Sehen Sie dazu auch

Überblick über die Pumpenkopfteile (> Seite 194)

7.5.1 Überblick über die Pumpenkopfteile



Die Abbildung zeigt die wichtigsten Teile des Pumpenkopfs.

Abbildung 89: Überblick über die Pumpenkopfteile

Nr.	Beschreibung
1	Pumpenkopf
2	Einlasskugelventil
3	Kolbendichtungen
4	Stützringe
5	Pumpenkopfdichtungen, für Schrauben der Hinterspülplatte (O-Ring- Dichtungen 6,5x1)
6	Pumpenkopfdichtungen, für Hinterspülplatten (O-Ring-Dichtungen 11x1,5)
7	Hinterspülplatten
8	Schrauben, Hinterspülplatten
9	Hinterspüldichtungen
10	Kolben
11	Schlauchverbinder für Hinterspülschläuche (schraubbar)
12	Auslasskugelventil
13	U-Rohr

7.5.2 Tauschen des Pumpenkopfs

Wann

Beschädigung des Pumpenkopfs

Gehen Sie wie folgt vor

- Entfernen Sie den Pumpenkopf. Siehe Entfernen des Pumpenkopfs (▶ Seite 195).
- Bauen Sie den Pumpenkopf ein. Siehe Einbauen des Pumpenkopfs (> Seite 197).

7.5.2.1 Entfernen des Pumpenkopfs

Wann

- Zum Tausch des Pumpenkopfs
- Bei der Wartung von Pumpenkopfteilen

Erforderliche Werkzeuge und zusätzliche Utensilien

- Sechskantschlüssel, Größe 6
- Lösungsmittel, geeignet zum Purgen, um schädliche Substanzen herauszuspülen

Vorbereitungen

1. Spülen Sie schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.

Betroffene Teile am Pumpenkopf



Abbildung 90: Pumpenkopf

Nr.	Beschreibung
1	Drucksensorkabel
2	Lösungsmittelschlauch am Einlasskugelventil
3	Pumpenkopfschraube

Nr.	Beschreibung
4-6	Hinterspülschläuche
7	Verbindungskapillare vom Pumpenkopf zum Flowmeter

Gehen Sie wie folgt vor

- Koppeln Sie die Kolben für den Pumpenkopf ab, für den Sie die Schritte durchführen möchten, entweder über die Tastatur (Tasten SELECT und DOCK) oder über das ePanel im Chromatographie-Datensystem.
- Warten Sie, bis der Abkoppelvorgang beendet ist, ehe Sie fortfahren. Das Abkoppeln ist beendet, wenn die Pumpe gestoppt hat und die LED neben der Taste FLOW für den gewählten Pumpenkopf grün blinkt.
- 3. Stecken Sie das Drucksensorkabel am Port **Pressure** ab.
- 4. Entfernen Sie den Lösungsmittelschlauch am Einlasskugelventil.
- 5. Entfernen Sie Kapillare, die den Pumpenkopf mit dem Flowmeter verbindet.
- 6. Stecken Sie die Hinterspülschläuche an den Schlauchverbindern am Pumpenkopf ab. Entfernen Sie nicht die Schlauchverbinder.
- 7. Lösen Sie die Pumpenkopfschraube mit dem Sechskantschlüssel.
- 8. Halten Sie den Pumpenkopf mit einer Hand und entfernen Sie die Pumpenkopfschraube.
- 9. Entfernen Sie den Pumpenkopf, indem Sie ihn circa 10 mm (0,5") nach vorn ziehen, weg von der Frontblende, und nach unten kippen. Die Kolben werden normalerweise zusammen mit dem Pumpenkopf aus der Pumpe entfernt. Wenn ein Kolben in der Pumpe bleibt, kippen Sie ihn und ziehen Sie ihn mit den Fingern heraus.

Sehen Sie dazu auch

- Einbauen des Pumpenkopfs (Seite 197)
- Purgen (Übersicht) (▶ Seite 65)

7.5.2.2 Einbauen des Pumpenkopfs

Wann

- Beim Einbau eines Austausch-Pumpenkopfs
- Bei der Wartung von Pumpenkopfteilen

Erforderliche Teile

Pumpenkopf

Erforderliche Werkzeuge und zusätzliche Utensilien

- Sechskantschlüssel, Größe 6
- Lösungsmittel, geeignet für Ihre Anwendung

Betroffene Teile am Pumpenkopf



Abbildung 91: Pumpenkopf

Nr.	Beschreibung
1	Drucksensorkabel
2	Lösungsmittelschlauch am Einlasskugelventil
3	Pumpenkopfschraube
4-6	Hinterspülschläuche
7	Verbindungskapillare vom Pumpenkopf zum Flowmeter

Beachten Sie Folgendes

Situation	Beschreibung
Wiedereinbau eines Pumpenkopfs	Vertauschen Sie die Pumpenköpfe nicht. Beachten Sie die Kennzeichnung am Drucksensorkabel (A für den linken Pumpenkopf, B für den rechten Pumpenkopf).
	Wenn Sie einen Kolben wieder einsetzen müssen, siehe Einbauen der Kolben (> Seite 201).
Einbau eines neuen Pumpenkopfs (Austausch- Pumpenkopf)	Neue Pumpenköpfe können beim erstmaligen Einbau in einer beliebigen Position eingebaut werden. Sie haben keine Kennzeichnung am Drucksensorkabel. Zur leichten Identifizierung bei künftigen Wartungsarbeiten sollten Sie sie gegebenenfalls entsprechend ihrer Einbauposition kennzeichnen.

Gehen Sie wie folgt vor

- 1. Überprüfen Sie, dass die Kolben soweit wie möglich in den Pumpenkopf eingesetzt sind.
- 2. Setzen Sie den Pumpenkopf in die Pumpe ein und ziehen Sie die Pumpenkopfschraube mit dem Sechskantschlüssel fest.
- 3. Verbinden Sie die Hinterspülschläuche mit den Schlauchverbindern auf dem Pumpenkopf.
- 4. Schließen Sie die Kapillare an, die den Pumpenkopf mit dem Flowmeter verbindet.
- 5. Verbinden Sie den Lösungsmittelschlauch mit dem Einlasskugelventil.
- 6. Stecken Sie das Kabel des Drucksensors am Port Pressure an.
- Koppeln Sie die Kolben entweder über die Tastatur an (Tasten SELECT und DOCK) oder über das ePanel im Chromatographie-Datensystem.
- Warten Sie, bis der Ankoppelvorgang beendet ist, ehe Sie fortfahren. Das Ankoppeln ist beendet, wenn die Pumpe gestoppt hat und die LED neben der Taste FLOW für den gewählten Pumpenkopf aufhört, grün zu blinken.
- 9. Schalten Sie die Pumpe über den Hauptnetzschalter aus und wieder ein.
- 10. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.
- 11. Prüfen Sie alle Flussverbindungen am Pumpenkopf optisch auf Anzeichen von Undichtigkeit. Ziehen Sie undichte Verbindungen fest, wenn das System druckfrei ist.

12. Testen Sie die Hinterspülung auf Dichtigkeit. Siehe Testen der Hinterspülung auf Undichtigkeit (▶ Seite 279).

13. Abhängig von der Situation:

Situation	Beschreibung
Wiedereinbau eines Pumpenkopfs	Testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: E04 - Detaillierter Dichtigkeitstest (Pumpe) .
Einbau eines neuen Pumpenkopfs (Austausch-Pumpenkopf)	 Führen Sie einen Abgleich der Drucksensoren durch. Führen Sie das Skript aus: C05 - System-Drucksensoren abgleichen.
	 Testen Sie die Pumpe auf Dichtigkeit. F ühren Sie das Skript aus: E04 - Detaillierter Dichtigkeitstest (Pumpe).

Sehen Sie dazu auch

- Entfernen des Pumpenkopfs (Seite 195)
- Purgen (Übersicht) (> Seite 65)

7.5.3 Tauschen der Kolben

Wann

Beschädigung des Kolbens

Gehen Sie wie folgt vor

- 1. Entfernen Sie den Kolben. Siehe Entfernen der Kolben (> Seite 200).
- 2. Bauen Sie den Kolben ein. Siehe Einbauen der Kolben (> Seite 201).

7.5.3.1 Entfernen der Kolben

Wann

- Zum Tausch eines Kolbens
- Bei der Wartung von Pumpenkopfteilen

Vorbereitungen

Entfernen Sie den Pumpenkopf. Siehe Entfernen des Pumpenkopfs (
 Seite 195).

Gehen Sie wie folgt vor

Ziehen Sie den Kolben, den Sie entfernen möchten, mit den Fingern aus dem Pumpenkopf heraus. Wenn Sie den Kolben bei der Wartung von Pumpenkopfteilen entfernen, fassen Sie ihn am Kolbenfuß an und vermeiden Sie es, die Kolbenoberfläche selbst zu berühren.



Abbildung 92: Entfernen der Kolben

Sehen Sie dazu auch

Einbauen der Kolben (▶ Seite 201)

7.5.3.2 Einbauen der Kolben

Wann

- Einbau eines Austauschkolbens, wenn Sie zuvor den Kolben entfernt hatten
- Bei der Wartung von Pumpenkopfteilen

Erforderliche Teile und zusätzliche Utensilien

- Kolben
- Isopropanol

Vorbereitungen

Wen Sie einen Kolben wiedereinbauen, den Sie entfernt haben, reinigen Sie den Kolben. Siehe Reinigen des Kolbens (▶ Seite 202).

Gehen Sie wie folgt vor



Abbildung 93: Einbauen der Kolben

Nr.	Beschreibung
1	Kolbenaufnahme
2	Kolbenfuß

- Pipettieren Sie einige Tropfen Isopropanol in die Kolbenaufnahme, in die Sie den Kolben einsetzen möchten.
 Wenn Sie einen Kolben wiedereinbauen, setzen Sie ihn an der Position ein, von der Sie ihn entfernt hatten.
- 2. Halten Sie den Kolben am Kolbenfuß.
- 3. Schieben Sie den Kolben soweit wie es geht in die Kolbenaufnahme hinein.
- Bauen Sie den Pumpenkopf ein. Siehe Einbauen des Pumpenkopfs (▶ Seite 197).

Sehen Sie dazu auch

Entfernen der Kolben (> Seite 200)

7.5.4 Reinigen des Kolbens

Wann

Bevor Sie einen Kolben wiedereinbauen, den Sie ausgebaut hatten

Erforderliche Utensilien

- Trockenes, fusselfreies Tuch
- Isopropanol

Gehen Sie wie folgt vor

- 1. Prüfen Sie den Kolben optisch auf Anzeichen einer Beschädigung.
- 2. Spülen Sie den Kolben vorsichtig mit Isopropanol.
- 3. Reiben Sie den Kolben mehrfach mit einem trocknen, fusselfreien Tuch ab.
- Halten Sie den Kolben fest und fahren Sie mit dem Fingernagel vorsichtig über die Kolbenoberfläche. Sie sollten keine unregelmäßigen oder rauen Stellen spüren.
- 5. Abhängig von der Situation:

Situation	Vorgehen
Kein sichtbarer Schaden Keine unregelmäßigen oder rauen Stellen	1. Spülen Sie den Kolben nochmals vorsichtig mit Isopropanol.
	 Reiben Sie den Kolben mehrfach mit einem trocknen, fusselfreien Tuch ab. Bauen Sie den Kolben wieder ein.
Kolben ist beschädigt oder verkratzt	Bauen Sie einen neuen Kolben ein.

Sehen Sie dazu auch

Einbauen der Kolben (> Seite 201)

7.5.5 Tauschen des Stützrings

Wann

Beschädigung des Stützrings

Gehen Sie wie folgt vor

- Entfernen Sie den Stützring. Siehe Entfernen des Stützrings (▶ Seite 203).
- Bauen Sie den Stützring ein. Siehe Einbauen des Stützrings (▶ Seite 205).

7.5.5.1 Entfernen des Stützrings

Wann

Zum Tausch des Stützrings

Erforderliche Werkzeuge

- Schraubendreher, Torx T30
- Dichtringwerkzeug



Abbildung 94: Dichtringwerkzeug (links: Werkzeug geschlossen; rechts: Werkzeug offen)

Nr.	Beschreibung
1	Ausbaupin
2	Einbaupin
3	Rändel zum Öffnen und Schließen des Dichtringwerkzeugs (Öffnen und Schließen ist nur bei Verwendung des Ausbaupins erforderlich)

Vorbereitungen

- Entfernen Sie den Pumpenkopf. Siehe Entfernen des Pumpenkopfs (> Seite 195).
- 2. Entfernen Sie den Kolben, für den Sie den Stützring ersetzen möchten. Siehe Entfernen der Kolben (▶ Seite 200).

Betroffene Teile am Pumpenkopf



Abbildung 95: Stützring (hier: für linken Kolben)

Nr.	Beschreibung
1	Pumpenkopf
2	Stützring
3	Pumpenkopfdichtungen (für Hinterspülplatte und Schrauben der Hinterspülplatte)
4	Hinterspülplatte
5	Schrauben, Hinterspülplatte (4 Schrauben)
6	Kolben

Gehen Sie wie folgt vor

1. Entfernen Sie mit dem Schraubendreher die Schrauben der Hinterspülplatte.



Abbildung 96: Schrauben der Hinterspülplatte (hier: linke Platte)

- 2. Nehmen Sie die Hinterspülplatte vom Pumpenkopf ab.
- Sollten sich eine oder mehrere Pumpenkopfdichtungen lösen, wenn Sie die Hinterspülplatte entfernen, heben Sie diese zum Wiedereinbau auf.

4. Entfernen Sie den Stützring: Drücken Sie den *Einbau*pin des Dichtringwerkzeugs leicht schräg in den Pumpenkopf. Verwenden Sie nicht den *Ausbau*pin.



Abbildung 97: Stützring (hier: für linken Kolben)

5. Entfernen Sie das Werkzeug. Der Stützring wird zusammen mit dem Werkzeug entfernt.

Sehen Sie dazu auch

Einbauen des Stützrings (Seite 205)

7.5.5.2 Einbauen des Stützrings

Wann

Einbau eines Austauschstützrings, wenn Sie zuvor den Stützring entfernt hatten

Erforderliche Teile

- Stützring
- Pumpenkopfdichtungen zum Wiedereinbau, falls erforderlich

Erforderliche Werkzeuge

- Schraubendreher, Torx T30
- Dichtringwerkzeug



Abbildung 98: Dichtringwerkzeug (links: Werkzeug geschlossen; rechts: Werkzeug offen)

Nr.	Beschreibung
1	Ausbaupin
2	Einbaupin
3	Rändel zum Öffnen und Schließen des Dichtringwerkzeugs (Öffnen und Schließen ist nur bei Verwendung des Ausbaupins erforderlich)

Betroffene Teile am Pumpenkopf



Abbildung 99: Stützring (hier: für linken Kolben)

Nr.	Beschreibung
1	Pumpenkopf
2	Stützring
3	Pumpenkopfdichtungen (für Hinterspülplatte und Schrauben der Hinterspülplatte)
4	Hinterspülplatte
5	Schrauben, Hinterspülplatte (4 Schrauben)
6	Kolben

Gehen Sie wie folgt vor

1. Schieben Sie den Stützring auf den *Einbau*pin des Dichtringwerkzeugs.



Abbildung 100: Aufschieben des Stützrings auf den Einbaupin

Nr.	Beschreibung
1	Stützring
2	Einbaupin, Dichtringwerkzeug

2. Drücken Sie das Dichtringwerkzeug mit dem Einbaupin in den Pumpenkopf und entfernen Sie das Werkzeug. Der Stützring bleibt im Pumpenkopf.



Abbildung 101: Stützring eingebaut (hier: für linken Kolben)

- Überprüfen Sie, ob die Pumpenkopfdichtungen korrekt eingesetzt sind. Wenn sich die Dichtungen gelöst haben oder nicht richtig in der Aufnahmenut liegen, setzen Sie die Dichtungen erneut ein. Siehe Einbauen der Pumpenkopfdichtungen (▶ Seite 221).
- 4. Legen Sie die Hinterspülplatte auf den Pumpenkopf und setzen Sie die Schrauben ein.



Abbildung 102: Schrauben der Hinterspülplatte (hier: linke Platte)

- 5. Ziehen Sie die Schrauben über Kreuz mit dem Schraubendreher fest.
- 6. Ziehen Sie die Schrauben nochmals über Kreuz fest.
- 7. Bauen Sie den Kolben ein. Siehe Einbauen der Kolben (> Seite 201).
- Bauen Sie den Pumpenkopf ein. Siehe Einbauen des Pumpenkopfs (▶ Seite 197).

Sehen Sie dazu auch

Entfernen des Stützrings (Seite 203)

7.5.6 Tauschen der Kolbendichtung

Wann

- Beschädigung der Kolbendichtung, undichte Dichtung
- Instabile Flussrate oder Basislinienrauschen

Gehen Sie wie folgt vor

- Entfernen Sie die Kolbendichtung. Siehe Entfernen der Kolbendichtung (> Seite 208).
- Bauen Sie die Kolbendichtung ein. Siehe Einbauen der Kolbendichtung (▶ Seite 211).

7.5.6.1 Entfernen der Kolbendichtung

Wann

Zum Tausch der Kolbendichtung

Erforderliche Werkzeuge

- Schraubendreher, Torx T30
- Dichtringwerkzeug



Abbildung 103: Dichtringwerkzeug (links: Werkzeug geschlossen; rechts: Werkzeug offen)

Nr.	Beschreibung
1	Ausbaupin
2	Einbaupin
3	Rändel zum Öffnen und Schließen des Dichtringwerkzeugs (Öffnen und Schließen ist nur bei Verwendung des Ausbaupins erforderlich)

Vorbereitungen

- Entfernen Sie den Pumpenkopf. Siehe Entfernen des Pumpenkopfs (▶ Seite 195).
- 2. Entfernen Sie den Kolben, für den Sie die Kolbendichtung ersetzen möchten. Siehe Entfernen der Kolben (▶ Seite 200).

Betroffene Teile am Pumpenkopf



Abbildung 104: Kolbendichtung (hier: für linken Kolben)

Nr.	Beschreibung
1	Pumpenkopf
2	Kolbendichtung
3	Stützring
4	Pumpenkopfdichtungen (für Hinterspülplatte und Schrauben der Hinterspülplatte)
5	Hinterspülplatte
6	Schrauben, Hinterspülplatte (4 Schrauben)
7	Kolben

Gehen Sie wie folgt vor

1. Entfernen Sie mit dem Schraubendreher die Schrauben der Hinterspülplatte.



Abbildung 105: Schrauben der Hinterspülplatte (hier: linke Platte)

- 2. Nehmen Sie die Hinterspülplatte vom Pumpenkopf ab.
- Sollten sich eine oder mehrere Pumpenkopfdichtungen lösen, wenn Sie die Hinterspülplatte entfernen, heben Sie diese zum Wiedereinbau auf.

4. Entfernen Sie den Stützring: Drücken Sie den *Einbau*pin des Dichtringwerkzeugs leicht schräg in den Pumpenkopf. Verwenden Sie nicht den *Ausbau*pin.



Abbildung 106: Stützring (hier: für linken Kolben)

- Entfernen Sie das Werkzeug. Der Stützring wird zusammen mit dem Werkzeug entfernt.
 Heben Sie den Stützring zum Wiedereinbau auf. Der Stützring ist kein Verschleißteil und muss nicht bei jedem Tausch der Kolbendichtung ersetzt werden.
- 6. Öffnen Sie das Dichtringwerkzeug und drücken Sie den *Ausbau*pin bis zum Anschlag in den Pumpenkopf.



Abbildung 107: Kolbendichtung (hier: für linken Kolben)

- Schließen Sie das Dichtringwerkzeug und entfernen Sie das Werkzeug aus dem Pumpenkopf. Die Dichtung wird zusammen mit dem Werkzeug entfernt.
- 8. Zum Entfernen der Dichtung vom Dichtringwerkzeug: Öffnen Sie das Werkzeug und ziehen Sie die Dichtung vom Werkzeug herunter. Eine Dichtung, die entfernt wurde, kann nicht wiederverwendet werden.

Sehen Sie dazu auch

Einbauen der Kolbendichtung (> Seite 211)

7.5.6.2 Einbauen der Kolbendichtung

Wann

Einbau einer Austausch-Kolbendichtung, wenn Sie die Dichtung zuvor entfernt hatten

Erforderliche Teile

- Kolbendichtung
- Stützring, der beim Entfernen der Kolbendichtung entfernt wurde Der Stützring ist kein Verschleißteil und kann wiedereingebaut werden. Bauen Sie nur dann einen neuen Stützring ein, wenn der Ring beschädigt ist.
- Pumpenkopfdichtungen zum Wiedereinbau, falls erforderlich

Erforderliche Werkzeuge und zusätzliche Utensilien

- Schraubendreher, Torx T30
- Spritze, gefüllt mit Isopropanol
- Reinigungsstäbchen
- Fusselfreies Tuch
- Dichtringwerkzeug



Abbildung 108: Dichtringwerkzeug (links: Werkzeug geschlossen; rechts: Werkzeug offen)

Nr.	Beschreibung
1	Ausbaupin
2	Einbaupin
3	Rändel zum Öffnen und Schließen des Dichtringwerkzeugs (Öffnen und Schließen ist nur bei Verwendung des Ausbaupins erforderlich)

Betroffene Teile am Pumpenkopf



Abbildung 109: Kolbendichtung (hier: für linken Kolben)

Nr.	Beschreibung
1	Pumpenkopf
2	Kolbendichtung
3	Stützring
4	Pumpenkopfdichtungen (für Hinterspülplatte und Schrauben der Hinterspülplatte)
5	Hinterspülplatte
6	Schrauben, Hinterspülplatte (4 Schrauben)
7	Kolben

Gehen Sie wie folgt vor

- 1. Reinigen Sie den Stützring von innen und außen mit Isopropanol, einem Reinigungsstäbchen und einem fusselfreien Tuch.
- 2. Pipettieren Sie einige Tropfen Isopropanol in die Aufnahme im Pumpenkopf, auf deren Rand die Kolbendichtung aufliegen wird.



Abbildung 110: Rand, auf dem Kolbendichtung aufliegt

3. Schieben Sie den Stützring und die Kolbendichtung auf den *Einbau*pin des Dichtringwerkzeugs. Die offene Seite der Kolbendichtung zeigt vom Werkzeug weg.



Abbildung 111: Aufschieben des Stützrings und der Kolbendichtung auf den Einbaupin

Nr.	Beschreibung
1	Kolbendichtung (offene Seite zeigt weg vom Werkzeug)
2	Stützring
3	Einbaupin, Dichtringwerkzeug

4. Drücken Sie das Dichtringwerkzeug mit dem Einbaupin bis zum Anschlag in den Pumpenkopf und entfernen Sie das Werkzeug. Die Kolbendichtung und der Stützring bleiben im Pumpenkopf.



Abbildung 112: Stützring eingebaut, verdeckt die Kolbendichtung (hier: für linken Kolben)

 Überprüfen Sie, ob die Pumpenkopfdichtungen korrekt eingesetzt sind. Wenn sich die Dichtungen gelöst haben oder nicht richtig in der Aufnahmenut liegen, setzen Sie die Dichtungen erneut ein. Siehe Einbauen der Pumpenkopfdichtungen (▶ Seite 221). 6. Legen Sie die Hinterspülplatte auf den Pumpenkopf und setzen Sie die Schrauben ein.



Abbildung 113: Schrauben der Hinterspülplatte (hier: linke Platte)

- 7. Ziehen Sie die Schrauben über Kreuz mit dem Schraubendreher fest.
- 8. Ziehen Sie die Schrauben nochmals über Kreuz fest.
- 9. Bauen Sie den Kolben ein. Siehe Einbauen der Kolben (> Seite 201).
- Bauen Sie den Pumpenkopf ein. Siehe Einbauen des Pumpenkopfs
 (▶ Seite 197). Bevor Sie jedoch das Skript zum Testen der Pumpe auf Dichtigkeit ausführen, betreiben Sie die Pumpe 30 Minuten lang bei einem Rückdruck von 80 MPa (empfohlen), damit die neuen Kolbendichtungen einlaufen können.

Sehen Sie dazu auch

Entfernen der Kolbendichtung (
 Seite 208)

7.5.7 Tauschen der Hinterspüldichtung

Wann

Beschädigung der Hinterspüldichtung, Flüssigkeit tritt unterhalb des Pumpenkopfs aus (undichte Dichtung)

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Hinterspüldichtung. Siehe Entfernen der Hinterspüldichtung (▶ Seite 215).
- Bauen Sie die Hinterspüldichtung ein. Siehe Einbauen der Hinterspüldichtung (▶ Seite 217).

7.5.7.1 Entfernen der Hinterspüldichtung

Wann

Zum Tausch der Hinterspüldichtung

Erforderliche Werkzeuge

Dichtringwerkzeug



Abbildung 114: Dichtringwerkzeug (links: Werkzeug geschlossen; rechts: Werkzeug offen)

Nr.	Beschreibung
1	Ausbaupin
2	Einbaupin
3	Rändel zum Öffnen und Schließen des Dichtringwerkzeugs (Öffnen und Schließen ist nur bei Verwendung des Ausbaupins erforderlich)

Vorbereitungen

- Entfernen Sie den Pumpenkopf. Siehe Entfernen des Pumpenkopfs (> Seite 195).
- 2. Entfernen Sie den Kolben, für den Sie die Hinterspüldichtung ersetzen möchten. Siehe Entfernen der Kolben (> Seite 200).

Betroffene Teile am Pumpenkopf



Abbildung 115: Hinterspüldichtung (hier: für linken Kolben)

Nr.	Beschreibung
1	Pumpenkopf
2	Hinterspülplatte
3	Hinterspüldichtung
4	Kolben

Gehen Sie wie folgt vor

1. Öffnen Sie das Dichtringwerkzeug und drücken Sie den *Ausbau*pin bis zum Anschlag in die Hinterspülplatte.



Abbildung 116: Hinterspüldichtung (hier: für linken Kolben)

- Schließen Sie das Dichtringwerkzeug und entfernen Sie das Werkzeug aus der Hinterspülplatte. Die Dichtung wird zusammen mit dem Werkzeug entfernt.
- 3. Zum Entfernen der Dichtung vom Dichtringwerkzeug: Öffnen Sie das Werkzeug und ziehen Sie die Dichtung vom Werkzeug herunter. Eine Dichtung, die entfernt wurde, kann nicht wiederverwendet werden.

Sehen Sie dazu auch

Einbauen der Hinterspüldichtung (> Seite 217)
7.5.7.2 Einbauen der Hinterspüldichtung

Wann

Einbau einer Austausch-Hinterspüldichtung, wenn Sie die Dichtung zuvor entfernt hatten

Erforderliche Teile und Werkzeuge

- Hinterspüldichtung
- Dichtringwerkzeug



Abbildung 117: Dichtringwerkzeug (links: Werkzeug geschlossen; rechts: Werkzeug offen)

Nr.	Beschreibung
1	Ausbaupin
2	Einbaupin
3	Rändel zum Öffnen und Schließen des Dichtringwerkzeugs (Öffnen und Schließen ist nur bei Verwendung des Ausbaupins erforderlich)

Betroffene Teile am Pumpenkopf





Nr.	Beschreibung
1	Pumpenkopf
2	Hinterspülplatte
3	Hinterspüldichtung
4	Kolben

 Schieben Sie die Hinterspüldichtung auf den Einbaupin des Dichtringwerkzeugs. Beachten Sie die Ausrichtung des Dichtrings.



Abbildung 119: Aufschieben der Hinterspüldichtung auf den Einbaupin

Nr.	Beschreibung
1	Hinterspüldichtung (der Pfeil zeigt an, welche Seite zum Werkzeug zeigt)
2	Einbaupin, Dichtringwerkzeug

2. Drücken Sie das Dichtringwerkzeug mit dem Einbaupin in die Hinterspülplatte und entfernen Sie das Werkzeug. Die Hinterspüldichtung bleibt in der Hinterspülplatte.



Abbildung 120: Hinterspüldichtung eingebaut (hier: für linken Kolben)

- 3. Bauen Sie den Kolben ein. Siehe Einbauen der Kolben (> Seite 201).
- Bauen Sie den Pumpenkopf ein. Siehe Einbauen des Pumpenkopfs (▶ Seite 197).

Sehen Sie dazu auch

Entfernen der Hinterspüldichtung (> Seite 215)

7.5.8 Tauschen der Pumpenkopfdichtungen

Wann

- Beschädigung der Pumpenkopfdichtungen
- Flüssigkeit tritt unterhalb des Pumpenkopfs aus (undichte Dichtung)

Gehen Sie wie folgt vor

- Entfernen Sie die Pumpenkopfdichtungen. Siehe Entfernen der Pumpenkopfdichtungen (▶ Seite 219).
- Bauen Sie die Pumpenkopfdichtungen ein. Siehe Einbauen der Pumpenkopfdichtungen (▶ Seite 221).

7.5.8.1 Entfernen der Pumpenkopfdichtungen

Wann

Zum Tausch der Pumpenkopfdichtungen

Erforderliche Werkzeuge

- Schraubendreher, Torx T30
- Pinzette

Vorbereitungen

- Entfernen Sie den Pumpenkopf. Siehe Entfernen des Pumpenkopfs (▶ Seite 195).
- 2. Entfernen Sie die Kolben. Siehe Entfernen der Kolben (> Seite 200).

Betroffene Teile am Pumpenkopf



Abbildung 121: Pumpenkopfdichtungen (hier: linke Seite des Pumpenkopfs)

Nr.	Beschreibung
1	Pumpenkopf
2	Pumpenkopfdichtungen, für Schrauben der Hinterspülplatte (O-Ring- Dichtungen 6,5x1)
3	Pumpenkopfdichtung, für Hinterspülplatte (O-Ring-Dichtung 11x1,5)
4	Hinterspülplatte

Nr.	Beschreibung
5	Schrauben der Hinterspülplatte
6	Kolben

Gehen Sie wie folgt vor

Empfohlen: Tauschen Sie alle Pumpenkopfdichtungen des Pumpenkopfs zur gleichen Zeit (linke Seite und rechte Seite des Pumpenkopfs).

1. Entfernen Sie mit dem Schraubendreher die Schrauben der Hinterspülplatte.



Abbildung 122: Schrauben der Hinterspülplatte (hier: linke Platte)

- 2. Nehmen Sie die Hinterspülplatte vom Pumpenkopf ab.
- 3. Entfernen Sie alle Pumpenkopfdichtungen. Das geht einfacher, wenn Sie eine Pinzette benutzen.



Abbildung 123: Pumpenkopfdichtungen (hier: linke Seite des Pumpenkopfs)

Nr.	Beschreibung
1	Pumpenkopfdichtungen, für Schrauben der Hinterspülplatte (O-Ring- Dichtungen 6,5x1), befinden sich im Pumpenkopf
2	Pumpenkopfdichtung, für Hinterspülplatte (O-Ring-Dichtung 11x1,5), befindet sich in der Hinterspülplatte, auf der Seite, die zum Pumpenkopf zeigt

Sehen Sie dazu auch

Einbauen der Pumpenkopfdichtungen () Seite 221)

7.5.8.2 Einbauen der Pumpenkopfdichtungen

Wann

- Einbau der Austausch-Pumpenkopfdichtungen, wenn Sie die Dichtungen zuvor entfernt hatten
- Wiedereinbau einer Pumpenkopfdichtung, die sich bei der Wartung an Pumpenkopfteilen gelöst hatte

Erforderliche Teile und Werkzeuge

- Pumpenkopfdichtungen
- Schraubendreher, Torx T30
- Spritze, gefüllt mit Isopropanol
- Fusselfreies Tuch

Betroffene Teile am Pumpenkopf



Abbildung 124: Pumpenkopfdichtungen (hier: linke Seite des Pumpenkopfs)

Nr.	Beschreibung
1	Pumpenkopf
2	Pumpenkopfdichtungen, für Schrauben der Hinterspülplatte (O-Ring- Dichtungen 6,5x1)
3	Pumpenkopfdichtung, für Hinterspülplatte (O-Ring-Dichtung 11x1,5)
4	Hinterspülplatte
5	Schrauben der Hinterspülplatte
6	Kolben

Gehen Sie wie folgt vor

- 1. Nur beim Wiedereinbau einer Pumpenkopfdichtung: Reinigen Sie die Pumpenkopfdichtung mit Isopropanol und einem fusselfreien Tuch.
- 2. Legen Sie die Pumpenkopfdichtung auf die Aufnahmenut.



Abbildung 125: Aufnahmenuten für Pumpenkopfdichtungen (hier: linke Seite des Pumpenkopfs), im Pumpenkopf und in der Hinterspülplatte

Nr.	Beschreibung
1	Aufnahmenuten für Pumpenkopfdichtungen im Pumpenkopf, Dichtungen für Schrauben der Hinterspülplatte (O-Ring-Dichtungen 6,5x1)
2	Aufnahmenut für Pumpenkopfdichtung in Hinterspülplatte (Seite, die zum Pumpenkopf zeigt), Dichtung für Hinterspülplatte (O-Ring-Dichtung 11x1,5)

3. Drücken Sie die Pumpenkopfdichtung mit den Fingern so weit wie möglich in die Nut.



Abbildung 126: Pumpenkopfdichtungen eingebaut, in Pumpenkopf und Hinterspülplatte (hier: linke Seite des Pumpenkopfs) 4. Legen Sie die Hinterspülplatte auf den Pumpenkopf und setzen Sie die Schrauben ein.



Abbildung 127: Schrauben der Hinterspülplatte (hier: linke Platte)

- 5. Ziehen Sie die Schrauben über Kreuz mit dem Schraubendreher fest.
- 6. Ziehen Sie die Schrauben nochmals über Kreuz fest.
- 7. Bauen Sie den Kolben ein. Siehe Einbauen der Kolben (> Seite 201).
- Bauen Sie den Pumpenkopf ein. Siehe Einbauen des Pumpenkopfs (> Seite 197).

Sehen Sie dazu auch

Entfernen der Pumpenkopfdichtungen (> Seite 219)

7.5.9 Warten der Pumpenkopf-Kugelventile

Der Pumpenkopf hat zwei Kugelventile: ein Einlasskugelventil und ein Auslasskugelventil.



Abbildung 128: Pumpenkopf-Kugelventile

Nr.	Beschreibung
1	Einlasskugelventil
2	Auslasskugelventil

Folgen Sie den Schritten für die Wartungsprozedur, die Sie ausführen möchten.

- Um das Einlasskugelventil zu tauschen, siehe Tauschen des Einlasskugelventils (Pumpenkopf) (▶ Seite 224).
- Um das Auslasskugelventil zu tauschen, siehe Tauschen des Auslasskugelventils (Pumpenkopf) (> Seite 226).
- Um die Kugelventile zu reinigen, siehe Reinigen der Kugelventile (Pumpenkopf) (▶ Seite 227).

TIPP Für das Kugelventil steht auch ein Skript zur Verfügung, dass Sie durch die Wartung führt. Führen Sie das Skript aus: **C10 – Pumpenkopf-Kugelventile reinigen oder auswechseln**.

7.5.9.1 Tauschen des Einlasskugelventils (Pumpenkopf)

Wann

- Kugelventilproblem, das nicht durch Reinigen behoben werden konnte (zum Beispiel, anhaltende Undichtigkeit)
- Beschädigung des Einlasskugelventils

Erforderliche Teile

Einlasskugelventil

Erforderliche Werkzeuge und zusätzliche Utensilien

- Lösungsmittel, geeignet zum Purgen, um schädliche Substanzen herauszuspülen
- Lösungsmittel, geeignet für Ihre Anwendung

Betroffene Teile am Pumpenkopf



Abbildung 129: Einlasskugelventil am Pumpenkopf

Nr.	Beschreibung
1	Lösungsmittelschlauch
2	Einlasskugelventil
3	Pumpenkopf

Vorbereitungen

1. Spülen Sie schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.

Gehen Sie wie folgt vor

- 1. Entfernen Sie den Lösungsmittelschlauch am Einlasskugelventil.
- 2. Lösen und entfernen Sie das Einlasskugelventil mit den Fingern.
- 3. Drehen Sie das neue Einlasskugelventil mit den Fingern fest.
- 4. Schließen Sie den Lösungsmittelschlauch an das Einlasskugelventil an.
- 5. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.
- 6. Überprüfen Sie das Kugelventil und die Anschlüsse optisch auf Anzeichen von Undichtigkeit. Ziehen Sie undichte Verbindungen nur fest, wenn das System druckfrei ist.

7. Testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: **E04 - Detaillierter Dichtigkeitstest (Pumpe)**.

Sehen Sie dazu auch

Purgen (Übersicht) (> Seite 65)

7.5.9.2 Tauschen des Auslasskugelventils (Pumpenkopf)

Wann

- Kugelventilproblem, das nicht durch Reinigen behoben werden konnte (zum Beispiel, anhaltende Undichtigkeit)
- Beschädigung des Auslasskugelventils

Erforderliche Teile

Auslasskugelventil

Erforderliche Werkzeuge und zusätzliche Utensilien

- Lösungsmittel, geeignet zum Purgen, um schädliche Substanzen herauszuspülen
- Lösungsmittel, geeignet für Ihre Anwendung

Betroffene Teile am Pumpenkopf





Nr.	Beschreibung
1	Pumpenkopf
2	Auslasskugelventil
3	U-Rohr

Vorbereitungen

1. Spülen Sie schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.

Gehen Sie wie folgt vor

- 1. Entfernen Sie am Pumpenkopf das U-Rohr.
- 2. Lösen und entfernen Sie das Auslasskugelventil mit den Fingern.
- 3. Setzen Sie das neue Auslasskugelventil ein und drehen Sie es mit den Fingern fest.
- 4. Verbinden Sie das U-Rohr am Pumpenkopf.
- 5. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.
- Überprüfen Sie das Kugelventil und die Anschlüsse optisch auf Anzeichen von Undichtigkeit. Ziehen Sie undichte Verbindungen nur fest, wenn das System druckfrei ist.
- 7. Testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: **E04 - Detaillierter Dichtigkeitstest (Pumpe)**.

Sehen Sie dazu auch

Purgen (Übersicht) (> Seite 65)

7.5.9.3 Reinigen der Kugelventile (Pumpenkopf)

Die Reinigungsprozedur umfasst Folgendes:

- Ultraschallreinigung mit reinem Isopropanol
- Einlaufenlassen der Kugelventile in der Pumpe bei hohem Rückdruck

Wann

- Einmal im Monat
- Verschmutzte oder fehlerhafte Kugelventile

Erforderliche Utensilien

- Spritze (12 mL), gefüllt mit Isopropanol
- Kleines Becherglas mit Isopropanol
- Ultraschallbad
- Lösungsmittel, geeignet für Ihre Anwendung (zum Purgen)
- Lösungsmittel, geeignet für die Einlaufprozedur
 In den meisten Fällen können Sie die Lösungsmittel Ihrer Applikation verwenden. Damit es nicht zu elektrostatischer Aufladung in der Gegendruckkapillare kommt, sollte die 50/50-Mischung der Lösungs-

mittel A und B mindestens 20% Wasser enthalten. Bei wasserfreien Lösungsmitteln, wie Normal-Phase-Lösungsmitteln, verwenden Sie 80% Isopropanol in Wasser sowohl in Kanal A als auch in Kanal B.

 Geeignete Gegendruckkapillare, die circa 75 MPa bei einem Fluss von 50 μL/min mit Wasser erzeugt (zum Beispiel, nanoViper-Kapillare, Fused Silica, Innendurchmesser: 20 μm, Länge: 350 mm)

Vorbereitungen

Entfernen Sie die Kugelventile, die Sie reinigen möchten.

Empfohlen: Reinigen Sie die Einlasskugelventile und Auslasskugelventile der beiden Pumpenköpfe zur gleichen Zeit.

Gehen Sie wie folgt vor

- 1. Füllen Sie mit der Spritze Isopropanol in Flussrichtung in die Kugelventile.
- 2. Legen Sie die Kugelventile in das Becherglas. Reinigen Sie sie circa 10 Minuten im Ultraschallbad.
- 3. Spülen Sie die Kugelventile in Flussrichtung mit Isopropanol aus der Spritze.
- 4. Setzen Sie die Kugelventile ein und drehen Sie sie fest. Schließen Sie das U-Rohr oder den Lösungsmittelschlauch wieder an.
- 5. Installieren Sie die Gegendruckkapillare am Pumpenausgang. Führen Sie das offene Ende der Kapillare in den Abfallbehälter.
- Lassen Sie die Pumpe die geeigneten Lösungsmittel für die Einlaufprozedur fördern: Beginnen Sie mit 50% B und einer Flussrate von 10 μL/min. Erhöhen Sie die Flussrate, bis der Systemdruck bei circa 90 MPa liegt und lassen Sie die Pumpe 20 Minuten laufen.
- Prüfen Sie den Pumpenkopf und die Kugelventile optisch auf Anzeichen von Undichtigkeit. Ziehen Sie undichte Verbindungen nur fest, wenn das System druckfrei ist.
- 8. Entfernen Sie die Gegendruckkapillare und schließen Sie das System wieder entsprechend Ihrer Anwendung an.
- 9. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.
- 10. Testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: **E04 -**Detaillierter Dichtigkeitstest (Pumpe).

Sehen Sie dazu auch

- Tauschen des Einlasskugelventils (Pumpenkopf) (Seite 224)
- Tauschen des Auslasskugelventils (Pumpenkopf) (> Seite 226)

7.6 Wartung des Flowmeters

Dieser Abschnitt beschreibt, wie Teile des Flowmeters gewartet oder getauscht werden. Folgen Sie den Schritten für die Wartungsprozedur, die Sie ausführen möchten.

7.6.1 Tauschen des Flowmeter-Filters

Wann

Filter verstopft

Diagnose-Skript empfiehlt Filtertausch

Erforderliche Teile

Flowmeter-Filterkit

Erforderliche Werkzeuge und zusätzliche Utensilien

- Maulschlüssel, Größe 13 mm
- Lösungsmittel, geeignet zum Ausspülen schädlicher Substanzen
- Lösungsmittel, geeignet für Ihre Anwendung

Betroffene Teile





Nr.	Beschreibung
1	Pumpenkopf
2	Verbindungskapillare vom Pumpenkopf zum Flowmeter
3	Flowmeter-Filter

Vorbereitungen

- 1. Spülen Sie schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.
- 2. Bereiten Sie den Austauschfilter vor.

Vorbereiten des Austauschfilters



Abbildung 132: Zusammensetzen des Flowmeter-Filters

Nr.	Beschreibung
1	Filterschraube
2	Filterschraube, Dichtung
3	Filterfritte

- 1. Schieben Sie die Dichtung mit den Fingern auf die Filterschraube, bis die Dichtung auf der Aufnahmekante aufliegt.
- 2. Pipettieren Sie einen Tropfen Lösungsmittel in die Filterschraube und setzen Sie die Filterfritte ein. Das Lösungsmittel hält die Fritte in der Schraube.

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Kapillare, die den Pumpenkopf und das Flowmeter verbindet.
- 2. Lösen Sie die Filterschraube mit dem Maulschlüssel und entfernen Sie den Filter aus der Pumpe.
- 3. Setzen Sie den neuen Filter ein und ziehen Sie die Filterschraube mit dem Maulschlüssel fest.
- 4. Schließen Sie die Kapillare an, die den Pumpenkopf und das Flowmeter verbindet.

- 5. Purgen Sie die Pumpe und das Flowmeter mit dem Lösungsmittel Ihrer Anwendung.
- Prüfen Sie die Verbindungen auf Anzeichen von Undichtigkeit. Ziehen Sie undichte Verbindungen fest, wenn das System druckfrei ist.
- 7. Überprüfen Sie, dass der neue Filter korrekt funktioniert. Führen Sie das Skript aus: **D01 Systemrückdruck testen**.
- 8. Testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: **E04 - Detaillierter Dichtigkeitstest (Pumpe)**.

Sehen Sie dazu auch

Purgen (Übersicht) (▶ Seite 65)

7.6.2 Tauschen der Flowmeter-Abfallleitung

Wann

Beschädigung der Abfallleitung oder wenn diese blockiert ist

Erforderliche Teile und zusätzliche Utensilien

- Flowmeter-Abfallleitung
- Lösungsmittel, geeignet zum Ausspülen schädlicher Substanzen

Betroffene Teile



Abbildung 133: Flowmeter-Abfallleitung (links: Übersicht; rechts: Detailansicht)

Nr.	Beschreibung
1	Flowmeter-Abfallleitung
2	Detektoreinheit der Hinterspülung

Nr.	Beschreibung
3	Drainage-Ablauf
4	Einbauposition für die Detektoreinheit der Hinterspülung
5	Waste-Ablauf (in die Systemdrainage)
6	Nut in der Wand des Drainage-Ablaufs
7	Leaksensor

Vorbereitungen

- 1. Spülen Sie schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.
- 2. Für den leichteren Zugang zur Wand des Drainage-Ablaufs und zum Waste-Ablauf entfernen Sie die Detektoreinheit der Hinterspülung (Trichter und Detektor).

Entfernen der Detektoreinheit der Hinterspülung

Fassen Sie die Detektoreinheit Hinterspülung (Trichter und Detektor) am Trichter und ziehen Sie die Einheit nach oben. Der Detektor mit Schlauch und Kabel kann mit dem Trichter verbunden bleiben.





Nr.	Beschreibung
1	Trichter, Hinterspüldetektor
2	Hinterspüldetektor
3	Detektorkabel
4	Hinterspülschlauch

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Abfallleitung am Port Purge Out am Flowmeter und am Waste-Ablauf.
- 2. Schließen Sie die neue Abfallleitung am Port Purge Out an.
- Führen Sie die Leitung durch die Nut in der Wand des Drainage-Ablauf in den Waste-Ablauf.
 Um einen falschen Leakalarm zu vermeiden, muss die Abfallleitung richtig im Waste-Ablauf sitzen. Ist das nicht der Fall, könnte sich Flüssigkeit im Drainage-Ablauf sammeln und den Leaksensor aktivieren.
- 4. Setzen Sie die Detektoreinheit der Hinterspülung wieder ein.

Sehen Sie dazu auch

Purgen (Übersicht) (▶ Seite 65)

7.7 Wartung der Lösungsmittelschläuche

Dieser Abschnitt beschreibt, wie Lösungsmittelschläuche oder der Lösungsmittelfilter gewartet oder getauscht werden. Folgen Sie den Schritten für die Wartungsprozedur, die Sie ausführen möchten.

7.7.1 Entleeren der Lösungsmittelschläuche

Wann

Um die Lösungsmittelschläuche zu entleeren, zum Beispiel, für Wartungs- oder Serviceprozeduren

Erforderliche Utensilien

Lösungsmittel, geeignet zum Ausspülen schädlicher Substanzen falls erforderlich

Gehen Sie wie folgt vor

- 1. Spülen Sie gegebenenfalls schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.
- Schrauben Sie den Deckel des Lösungsmittelbehälters ab und entfernen Sie den Deckel zusammen mit dem Lösungsmittelschlauch vom Behälter.
- 3. Starten Sie einen Purge-Zyklus wie erforderlich.
- Stoppen Sie den Purge-Zyklus, sobald der Lösungsmittelschlauch leer ist.

Sehen Sie dazu auch

Purgen (Übersicht) (▶ Seite 65)

7.7.2 Tauschen der Lösungsmittelschläuche vom Behälter zur Pumpe

Wann

Beschädigung des Lösungsmittelschlauchs oder wenn dieser blockiert ist

Erforderliche Teile und zusätzliche Utensilien

- Lösungsmittelschläuche vom Behälter zur Pumpe
- Lösungsmittel, geeignet für Ihre Anwendung

Betroffene Teile



Abbildung 135: Lösungsmittelschläuche vom Behälter zur Pumpe

Nr.	Beschreibung
1	Lösungsmittelschläuche von den Behältern, zur Pumpe geführt durch alle Module oberhalb der Pumpe im Systemturm, mit
1a	Lösungsmittelschlauch, Kanal A
1b	Lösungsmittelschlauch, Kanal B
2	Absperrventile, A und B
3	Detektoreinheit der Hinterspülung

Vorbereitungen

- 1. Entleeren Sie die Lösungsmittelschläuche. Siehe Entleeren der Lösungsmittelschläuche (▶ Seite 234).
- Wenn Sie den Lösungsmittelfilter zusammen mit dem Lösungsmittelschlauch tauschen möchten, bereiten Sie den Filter vor. Siehe Tauschen des Lösungsmittelfilters (▶ Seite 240).
- 3. Entfernen Sie für den leichteren Zugang die Detektoreinheit der Hinterspülung (Trichter und Detektor).

Entfernen der Detektoreinheit der Hinterspülung

Fassen Sie die Detektoreinheit Hinterspülung (Trichter und Detektor) am Trichter und ziehen Sie die Einheit nach oben. Der Detektor mit Schlauch und Kabel kann mit dem Trichter verbunden bleiben.





Nr.	Beschreibung
1	Trichter, Hinterspüldetektor
2	Hinterspüldetektor
3	Detektorkabel
4	Hinterspülschlauch

Entfernen des Lösungsmittelschlauchs

- 1. Entfernen Sie den Lösungsmittelschlauch aus dem Behälterdeckel:
 - a) Entfernen Sie den Filter vom Lösungsmittelschlauch.
 - b) Entfernen Sie die Schlauchführung.
 - c) Ziehen Sie den Lösungsmittelschlauch aus dem Behälterdeckel heraus.
- 2. Lösen und entfernen Sie den Lösungsmittelschlauch in der Pumpe.
- Entfernen Sie den Lösungsmittelschlauch aus der Pumpe und aus allen Modulen, die sich oberhalb der Pumpe im Systemturm befinden. Wenn Sie den Lösungsmittelschlauch aus den Schlauchführungen entfernen, ziehen Sie nicht an den anderen Schläuchen in den Führungen.

Führen des neuen Lösungsmittelschlauchs zum Solvent Rack

- 1. Verbinden Sie jeden Lösungsmittelschlauch mit dem Eingangsport in der Pumpe, für den er gedacht ist (beachten Sie die Kennzeichnung an den Lösungsmittelschläuchen und in der Pumpe).
- Führen Sie die Lösungsmittelschläuche von der Pumpe zum Solvent Rack, durch die Schlauchführungen in der Pumpe und in allen Modulen, die sich oberhalb der Pumpe im Systemturm befinden. Beachten Sie die Hinweise zur Verlegung in Kapillar- und Schlauchführung durch das System (▶ Seite 119).
- 3. Führen Sie die Lösungsmittelschläuche durch das Führungsloch im Solvent Rack und drücken Sie sie in eine oder mehrere Schlauchführungen.





Nr.	Beschreibung
А	Führungsloch (Solvent Rack)
1-4	Schlauchführungen

Verbinden des Lösungsmittelschlauchs mit dem Lösungsmittelbehälter



Abbildung 138: Vorbereiten des Lösungsmittelschlauchs und des Behälterdeckels

Nr.	Beschreibung
1	Lösungsmittelschlauch
2	Schlauchführung

Nr.	Beschreibung
3	Behälterdeckel
4	Verschlusskappen
5	Lösungsmittelfilter mit Schlauchadapter

- 1. Führen Sie jeden Lösungsmittelschlauch durch eine Schlauchführung.
- 2. Führen Sie den Lösungsmittelschlauch durch eine Öffnung in einem Deckel für einen Lösungsmittelbehälter. Die Schlauchführung verhindert, dass der Lösungsmittelschlauch im Behälter verrutscht.
- 3. Verschließen Sie die Öffnungen in den Behälterdeckeln mit den Verschlusskappen.
- 4. Schieben Sie einen Lösungsmittelfilter auf jeden Lösungsmittelschlauch.
- 5. Drehen Sie den Flaschendeckel handfest an. Drücken Sie die Schlauchführung in die Öffnung im Behälterdeckel, um den Schlauch im Deckel zu arretieren.
- 6. Platzieren Sie die Lösungsmittelschläuche gerade in den Schlauchführungen.
- 7. Prüfen Sie die Lösungsmittelschläuche über den gesamten Flussweg hinweg. Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
- 8. Setzen Sie, falls erforderlich, die Detektoreinheit der Hinterspülung wieder ein.
- 9. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.

Sehen Sie dazu auch

Purgen (Übersicht) (▶ Seite 65)

7.7.3 Tauschen der Lösungsmittelschläuche vom Absperrventil zum Pumpenkopf

Wann

Beschädigung des Lösungsmittelschlauchs oder wenn dieser blockiert ist

Erforderliche Teile und zusätzliche Utensilien

- Lösungsmittelschläuche vom Absperrventil zum Pumpenkopf
- Lösungsmittel, geeignet zum Ausspülen schädlicher Substanzen
- Lösungsmittel, geeignet für Ihre Anwendung

Vorbereitungen

• Spülen Sie schädliche Substanzen aus, indem Sie die Pumpe mit einem geeigneten Lösungsmittel purgen.

Betroffene Teile



Abbildung 139: Lösungsmittelschläuche von den Absperrventilen zu den Pumpenköpfen

Nr.	Beschreibung
1	Absperrventile, A und B
2	Pumpenköpfe, A und B
3	Lösungsmittelschlauch vom Absperrventil B zum Pumpenkopf B
4	Lösungsmittelschlauch vom Absperrventil A zum Pumpenkopf A

Gehen Sie wie folgt vor

- 1. Entfernen Sie den Lösungsmittelschlauch am Absperrventil und am Pumpenkopf.
- 2. Schließen Sie den neuen Lösungsmittelschlauch an.
- 3. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.

Sehen Sie dazu auch

Purgen (Übersicht) (▶ Seite 65)

7.7.4 Tauschen des Lösungsmittelfilters

Wann

- Während regelmäßiger Wartungsarbeiten
- Durchlässigkeit der Fritte im Lösungsmittelfilter beeinträchtigt
- Lösungsmittelfilter beschädigt

Erforderliche Teile und zusätzliche Utensilien

- Wie zutreffend:
 - Filterfritte
 - Lösungsmittelfilter (mit Filterhalter, Filterfritte, und Schlauchadapter)
- Lösungsmittel, geeignet für Ihre Anwendung

Betroffene Teile



Abbildung 140: Lösungsmittelfilter (links: bei Tausch der Filterfritte; rechts: bei Tausch des Lösungsmittelfilters)

Nr.	Beschreibung
1	Schlauchadapter
2	Filterhalter, Oberteil
3	Filterfritte
4	Filterhalter, Unterteil

Vorbereitungen

- 1. Schrauben Sie den Deckel des Lösungsmittelbehälters ab und entfernen Sie den Deckel zusammen mit dem Lösungsmittelschlauch vom Behälter.
- 2. Entfernen Sie den Filterhalter (mit Adapter) vom Lösungsmittelschlauch.

Gehen Sie wie folgt vor

Folgen Sie den entsprechenden Schritten zum Tauschen der Filterfritte oder des Filters.

Schritte zum Tauschen der Fritte im Lösungsmittelfilter

- 1. Schrauben Sie dazu den Filterhalter auf und tauschen Sie die Fritte aus.
- 2. Setzen Sie den Lösungsmittelfilter zusammen; tragen Sie dabei geeignete saubere Handschuhe:
 - a) Legen Sie die Fritte waagrecht in den Filterhalter (Unterteil).
 - b) Schrauben Sie das Oberteil des Filters fest auf das Unterteil.
- 3. Schieben Sie den Filterhalter mit Adapter auf den Lösungsmittelschlauch.
- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 5. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung, um die Luft aus den Lösungsmittelschläuchen zu entfernen.

Schritte zum Tauschen des Lösungsmittelfilters

- 1. Setzen Sie den Lösungsmittelfilter zusammen; tragen Sie dabei geeignete saubere Handschuhe:
 - a) Setzen Sie den Schlauchadapter in den Filterhalter (Oberteil) ein.
 - b) Drehen Sie den Filterhalter (Oberteil) mit Adapter auf den Kopf und drücken Sie ihn gegen eine geeignete Oberfläche, so dass der Adapter fest im Filterhalter sitzt.
 - c) Legen Sie die Fritte waagrecht in den Filterhalter (Unterteil).
 - d) Schrauben Sie das Oberteil des Filters fest auf das Unterteil.
- 2. Schieben Sie den Filterhalter mit Adapter auf den Lösungsmittelschlauch.
- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 4. Purgen Sie die Pumpe mit dem Lösungsmittel Ihrer Anwendung.

Sehen Sie dazu auch

Purgen (Übersicht) (▶ Seite 65)

7.8 Wartung der Probenschleife

Wann

- Probenschleife ist verstopft oder beschädigt
- Anderes Probenschleifenvolumen benötigt

Erforderliche Teile

Ersatz-Probenschleife oder Probenschleife einer anderen Größe

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Probenschleife.
- 2. Bauen Sie die Ersatz-Probenschleife ein.
- Schließen Sie die Tauschprozedur ab und wählen Sie die entsprechende Probenschleife im Chromatographie-Datensystem aus.

TIPP Für die Probenschleife steht auch ein Skript zur Verfügung, dass Sie durch die Wartung führt. Wenn Sie Hilfe benötigen, führen Sie das Skript aus: **A07 - Probenschleife ändern**.

7.8.1 Komponenten der Probenschleife





Nr.	Beschreibung
1	Probenschleifen-Fitting an der Nadeleinheit
2	Probenschleife
3	Probenschleifen-Platte
4	Befestigungsschraube für die Probenschleifen-Platte
5	Aussparung in der Probenschleifen-Platte

7.8.2 Entfernen der Probenschleife

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Vorbereitungen

- 1. Fahren Sie die Nadeleinheit in die Service-Position.
- 2. Entfernen Sie die Isolierabdeckung.
- 3. Sichern Sie die Nadeleinheit.
- 4. Schalten Sie den Autosampler über den Hauptnetzschalter aus.

Gehen Sie wie folgt vor

- Lösen Sie mit dem Schraubendreher leicht die Befestigungsschraube. Die Schraube kann locker am Halter der Probenschleifeneinheit angeschraubt bleiben.
- 2. Entfernen Sie das Probenschleifen-Fitting von der Nadeleinheit. Verwenden Sie zum leichteren Entfernen den Drehmoment-Schraubendreher mit Viper-Steckaufsatz.
- 3. Entfernen Sie das Probenschleifen-Fitting vom Injektionsventil.
- Drücken Sie die Probenschleifeneinheit leicht nach oben, sodass die Befestigungsschraube in die große Öffnung der Probenschleifen-Platte passt. Entfernen Sie die Probenschleifeneinheit von der Nadeleinheit.



Abbildung 142: Entfernen der Probenschleife

 Wenn die Probenschleife wieder eingebaut werden soll: Schützen Sie offene Probenschleifen-Fittinge mit einer Schutzkappe.

7.8.3 Einbauen der Probenschleife

Erforderliche Teile und Werkzeuge

- *Wenn die Probenschleife getauscht werden soll:* Ersatz-Probenschleife
- Schraubendreher, Torx T10
- Drehmoment-Schraubendreher mit Verlängerung und Viper-Steckaufsatz

Gehen Sie wie folgt vor

- Entfernen Sie die Schutzkappe vom unteren Probenschleifen-Fitting und schließen Sie die Probenschleife an das Injektionsventil an. Weitere Informationen zu den Anschlüssen finden Sie unter Anschlussbelegung der Autosampler-Ventile (> Seite 48).
- 2. Entfernen Sie die Kappe vom oberen Probenschleifen-Fitting und schließen Sie die Probenschleife an die Nadeleinheit an.
- 3. Hängen Sie die Probenschleifen-Platte am Halter für die Probenschleifen-Einheit ein:
 - a) Hängen Sie die Schrauben-Aussparung in der Probenschleifen-Platte an die Befestigungsschraube. Achten Sie darauf, dass die Schraube gelockert ist.
 - b) Richten Sie die längliche Aussparung in der Platte an der Halterung aus.
- Führen Sie das Probenschleifen-Fitting in die Nadeleinheit und schrauben Sie es ein.
 Vergewissern Sie sich, dass die Kapillare von der Nadeleinheit wegführend so ausgerichtet ist, dass die Kapillare aus dem Stecker auf 5 Uhr zeigt.
- 5. Ziehen Sie das Probenschleifen-Fitting an der Nadeleinheit mit dem Drehmoment-Schraubendreher fest. Nach dem Festziehen sollte die Kapillare auf 6 Uhr zeigen.

- 6. Ziehen Sie die Befestigungsschraube für die Probenschleifen-Platte mit dem Schraubendreher fest.

Abbildung 143: Ausrichten der Probenschleifeneinheit an der Schraube

7. Führen Sie die Kapillare, die zum Injektionsventil führt, durch die Halterung der Isolierabdeckung.



Abbildung 144: Führen der Kapillare durch die Halterung der Isolierabdeckung

8. Schließen Sie die Tauschprozedur ab. Siehe Abschließen des Probenschleifen-Tauschs (▶ Seite 245).

7.8.4 Abschließen des Probenschleifen-Tauschs

Gehen Sie nach dem Einbau der Probenschleife wie folgt vor:

- 1. Entsichern Sie die Nadeleinheit.
- Prüfen Sie, dass sich die Probenschleife leichtgängig bewegt, indem Sie die Nadeleinheit in den Probenraum hinein schieben.
 Wenn die Probenschleife an der Innenwand festhängt, bauen Sie die Probenschleife erneut aus und ein entsprechend den Schritten in diesem Abschnitt. Achten Sie darauf, dass die Probenschleife korrekt ausgerichtet ist und der Stecker mit der Kapillare auf 6 Uhr zeigt.

- 3. Bauen Sie die Isolierabdeckung ein.
- 4. Schalten Sie den Autosampler über den Hauptnetzschalter ein. Der Autosampler führt einen Selbsttest durch. Die Nadel wird initialisiert und justiert ihre Position.
- 5. Falls erforderlich, stellen Sie den Parameter **Nominal Loop Volume** manuell für die eingebaute Probenschleife ein. Beachten Sie die Volumen, die auf dem Schild der Probenschleife angegeben sind.

Beschreibung	Nominalvolumen der Probenschleife	Gesamtvolumen der Probenschleife
Probenschleife, 10 μL	10 μL	23 μL
Probenschleife, 25 μL	25 μL	50 μL
Probenschleife, 100 μL	100 μL	130 μL

6. Bereiten Sie den Autosampler für den Betrieb vor. Siehe Vorbereiten des Systems für den Betrieb (▶ Seite 160).

7.9 Wartung der Nadeleinheit

7.9.1 Komponenten der Nadeleinheit und Nadelsitz



Abbildung 145: Komponenten der Nadeleinheit und Nadelsitz

Nr.	Beschreibung
1	Transportsicherungsschraube
2	Sperre für den Niederhalter
3	Positionierkante am Niederhalter der Nadeleinheit
4	Befestigungsschraube für den Niederhalter
5	Schrauben zur Befestigung der Nadeleinheit am Nadelschlitten
6	Probennadel
7	Niederhalter
8	Nadelsitz

7.9.2 Sichern der Nadeleinheit

Wann

Vor Wartungsmaßnahmen oder für den Transport

ACHTUNG

Wenn die Nadeleinheit während Wartungsarbeiten oder einem Transport nicht gesichert ist, kann sie den Autosampler beschädigen. Sichern Sie die Nadeleinheit, wenn angewiesen, um Schäden am Autosampler zu vermeiden.

Erforderliche Werkzeuge

Schlitzschraubendreher (optional)

Vorbereitungen

Entfernen Sie die Isolierabdeckung. Siehe Entfernen der Isolierabdeckung (▶ Seite 98).

Hochfahren der Nadeleinheit in Service-Position

Drücken Sie die Taste **SERVICE** auf der Tastatur einmal, um die Nadeleinheit aus dem Nadelsitz hoch in die erste Service-Position zu fahren.

Festziehen der Transportsicherungsschraube

Ziehen Sie die Sicherungsschraube oberhalb der Nadeleinheit mit den Fingern fest, bis sie sich nicht mehr weiter anziehen lässt. Die Verwendung eines Schlitzschraubendrehers erleichtert gegebenenfalls das Drehen der Schraube.



Abbildung 146: Festziehen der Transportsicherungsschraube

7.9.3 Entsichern der Nadeleinheit

Erforderliche Werkzeuge

Schlitzschraubendreher (optional)

Lösen der Transportsicherungsschraube

ACHTUNG

Wenn die Transportsicherungsschraube nicht korrekt gelöst wurde, funktioniert der Autosampler möglicherweise nicht richtig oder kann beschädigt werden. Vergewissern Sie sich immer, dass die Transportsicherungsschraube vollständig gelöst ist, bevor Sie den Autosampler einschalten.

Lösen Sie die Transportsicherungsschraube oberhalb der Nadeleinheit: Drehen Sie die unverlierbare Schraube gegen den Uhrzeigersinn mit Ihren Fingern, bis sie locker in ihrer Feder sitzt. Die Verwendung eines Schlitzschraubendrehers erleichtert gegebenenfalls das Drehen der Schraube.



Abbildung 147: Lösen der Transportsicherungsschraube

Prüfen der Nadeleinheit

Nach Erfordernis:

Zustand des Autosamplers	Schritte
Der Autosampler war während der Wartungsprozedur ausgeschaltet	Prüfen Sie, dass sich die Nadeleinheit leichtgängig bewegt: Schieben Sie die Nadeleinheit entlang des horizontalen Nadelantriebs in den Probenraum hinein. Siehe Bild unten. Wenn sich die Nadeleinheit leicht verschieben lässt, ist die Transportsicherungsschraube korrekt gelöst.
Der Autosampler ist während der Wartungsprozedur eingeschaltet geblieben	Drücken Sie die Taste SERVICE , um die Nadeleinheit in den Probenraum zu bewegen.



Abbildung 148: Verschieben der Nadeleinheit (bei ausgeschaltetem Autosampler)

Einrichten der Nadeleinheit nach der Service-Position

Nach Erfordernis:

Zustand des Autosamplers	Schritte
Der Autosampler war während der Wartungsprozedur ausgeschaltet	Die Nadeleinheit wird automatisch beim Einschalten des Autosamplers initialisiert.
Der Autosampler ist während der Wartungsprozedur eingeschaltet geblieben	Abhängig davon, welche Service-Position genutzt wurde:
	 Drücken Sie die Taste SERVICE und warten Sie, bis sich die Nadel in die erste Service-Position bewegt hat (Nadel ist oben).
	 Drücken Sie die Taste SERVICE erneut und warten Sie, bis sich die Nadel in die zweite Service-Position bewegt hat (Nadel über Waschport).
	 Drücken Sie die Taste SERVICE ein drittes Mal, um die Nadeleinheit nach der Service-Position zu initialisieren. Prüfen Sie, ob sich die Probenschleife frei bewegen kann.

7.9.4 Tauschen des Nadelsitzes und der Nadeleinheit

Wann

Je nach Art des Problems kann es erforderlich sein, Nadelsitz und Nadeleinheit zusammen oder getrennt zu tauschen.

Wann	Abhilfe
Nadelsitz ist verstopft	1. Führen Sie das Skript aus: C21 - Nadeleinheit und Nadelsitz reinigen oder tauschen
	 2. Wenn das Skript die Blockierung nicht entfernt: Tauschen Sie den Nadelsitz, siehe Tauschen des Nadelsitzes (▶ Seite 252).
	3. Überprüfen Sie die Nadeleinheit auf Schäden.
Nadelsitz ist undicht	1. Ziehen Sie den Nadelsitz mit dem Drehmoment- Schraubendreher mit Nadelsitz-Steckaufsatz fest. Siehe auch Prüfen der Autosampler-Ventile auf Dichtigkeit (▶ Seite 296).
	Wenn der Nadelsitz immer noch undicht ist: Tauschen Sie Nadelsitz und Nadeleinheit zusammen.
	2. Tauschen Sie den Nadelsitz, siehe Tauschen des Nadelsitzes () Seite 252).
	3. Tauschen Sie die Nadeleinheit, siehe Tauschen der Nadeleinheit (▶ Seite 255).
Nadel ist beschädigt	Tauschen Sie Nadelsitz und Nadeleinheit zusammen:
	1. Tauschen Sie den Nadelsitz, siehe Tauschen des Nadelsitzes () Seite 252).
	2. Tauschen Sie die Nadeleinheit, siehe Tauschen der Nadeleinheit (▶ Seite 255).

Erforderliche Teile

- Nadeleinheit
- Nadelsitz

Erforderliche Werkzeuge

- Schraubendreher, Torx T10
- Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz



Abbildung 149: Werkzeug, um den Nadelsitz anzuziehen

Nr.	Beschreibung
1	Nadelsitz-Steckaufsatz
2	Basis des Drehmoment-Schraubendrehers

TIPP Es ist möglich, ein Skript auszuführen, um Schritt-für-Schritt-Anweisungen direkt am Gerät zu erhalten. Für weitere Informationen führen Sie das Skript aus: **C21 - Nadeleinheit und Nadelsitz reinigen oder tauschen**.

7.9.4.1 Tauschen des Nadelsitzes

Gehen Sie wie folgt vor

- 1. Drücken Sie die Taste **SERVICE** zweimal, um die Nadel in die zweite Service-Position über dem Waschport zu bewegen.
- 2. Schalten Sie den Autosampler über den Hauptnetzschalter aus.
- 3. Entfernen Sie die Isolierabdeckung.
- 4. Entfernen Sie die Kapillare aus der Halterung der Isolierabdeckung.


5. Drücken Sie die Halterung der Isolierabdeckung zusammen und schieben Sie sie hoch, um sie zu entfernen.

Abbildung 150: Entfernen der Halterung der Isolierabdeckung

ACHTUNG—Der Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz kann beschädigt werden, wenn er nicht korrekt verwendet wird.

- Setzen Sie den Nadelsitz-Steckaufsatz vollständig in den Drehmoment-Schraubendreher ein.
- Setzen Sie den Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz direkt auf den Nadelsitz.
- Drücken oder ziehen Sie das Werkzeug nie mit unnötiger Kraft.
- 6. Setzen Sie den Nadelsitz-Steckaufsatz in den Drehmoment-Schraubendreher ein.
- 7. Setzen Sie den Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz auf den Nadelsitz.

8. Drehen Sie das Werkzeug gegen den Uhrzeigersinn, um den

Abbildung 151: Entfernen des Nadelsitzes mit dem Drehmoment-Schraubendreher

9. Entpacken Sie den Ersatz-Nadelsitz.

Nadelsitz zu entfernen.

10. Setzen Sie den Nadelsitz ein und drehen Sie ihm mit den Fingern im Uhrzeigersinn fest.



Abbildung 152: Einsetzen des Nadelsitzes

11. Verwenden Sie den Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz und drehen Sie ihn im Uhrzeigersinn, um den Nadelsitz festzuziehen, bis die Drehmomentbegrenzung auslöst (erkennbar am hörbaren Klicken).

ACHTUNG—Andere Werkzeuge können den Nadelsitz beschädigen

Der Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz ist zum Festziehen des Nadelsitzes gedacht. Bei Verwendung dieses Werkzeugs wird der Nadelsitz korrekt festgezogen und wird nicht zu fest angezogen. Verwenden Sie kein anderes Werkzeug.



Abbildung 153: Festziehen des Nadelsitzes

12. Je nach Art des Problems fahren Sie mit den Schritten in Kapitel Tauschen der Nadeleinheit (▷ Seite 255) oder Abschließen des Tauschs von Nadelsitz und Nadeleinheit (▷ Seite 258) fort.

7.9.4.2 Tauschen der Nadeleinheit

TIPP Eine beschädigte Nadeleinheit verursacht auch Schäden am Nadelsitz. Daher wird empfohlen, den Nadelsitz immer zusammen mit der Nadeleinheit zu tauschen.

Gehen Sie wie folgt vor

- Entfernen Sie die Probenschleife. Folgen Sie den Schritten zum Entfernen der Probenschleife in Kapitel Entfernen der Probenschleife (▶ Seite 243). Die Verbindung zum Injektionsventil kann angeschlossen bleiben.
- 2. Ziehen Sie die Nadeleinheit mit der Hand leicht nach vorn heraus.
- 3. Sichern Sie die Nadeleinheit.



4. Lösen Sie die zwei Schrauben, die die Nadeleinheit am Nadelschlitten befestigen, mit dem Schraubendreher.

Abbildung 154: Abschrauben der Nadeleinheit

Nr.	Beschreibung
1	Schrauben zur Befestigung der Nadeleinheit am Nadelschlitten

5. Ziehen Sie die Nadeleinheit leicht nach vorn und weg vom Nadelschlitten.

ACHTUNG—Die Öffnung für die Positionierkante ist empfindlich gegen Staub und Schmutzpartikel.

Achten Sie darauf, dass kein Staub und keine Schmutzpartikel in das Innere des Autosamplers gelangen.



WARNUNG—Scharfe Spitze der Injektionsnadel

Die Injektionsnadel besitzt eine scharfe Spitze, die Verletzungen der Haut verursachen kann.

Berühren Sie niemals die Nadelspitze, um Personenschäden zu vermeiden.

ACHTUNG—Die Nadel in der Nadeleinheit ist nicht befestigt.

- Entfernen Sie die Nadeleinheit vorsichtig, nachdem Sie sie abgeschraubt haben.
- Gehen Sie vorsichtig mit der Ersatz-Nadeleinheit um, insbesondere nach dem Entfernen der Transportsicherung.
- 6. Entpacken Sie die Ersatz-Nadeleinheit.
- Verwenden Sie die Nadeleinheit mit montierter Transportsicherung. Die Transportsicherung wird beim Einsetzen der Nadeleinheit automatisch gelöst und kann anschließend entfernt werden.
- 8. Positionieren Sie die Ersatz-Nadeleinheit auf dem Nadelschlitten. Beachten Sie folgende Hinweise:
 - Die vorstehende Kante auf der Rückseite der Nadeleinheit muss in die vorgesehene Öffnung am Nadelschlitten passen (siehe Abbildung).
 - Die Erhöhung an der Rückseite der Nadeleinheit muss in die Aussparung am Nadelschlitten passen.



Abbildung 155: Montieren der Nadeleinheit am Nadelschlitten

Nr.	Beschreibung
1	Öffnung am Nadelschlitten
2	Vorstehende Kante an der Nadeleinheit

9. Setzen Sie die zwei Schrauben an der Nadeleinheit ein und ziehen Sie sie fest.



10. Prüfen Sie, ob sich Sperre und Niederhalter leicht anheben lassen.

Abbildung 156: Anheben der Sperre für den Niederhalter

Nr.	Beschreibung
1	Sperre für den Niederhalter
2	Niederhalter

- 11. Entsichern Sie die Nadeleinheit.
- 12. Bewegen Sie die Nadeleinheit leicht nach hinten über den Waschport.
- 13. Fahren Sie mit den Schritten in Kapitel Abschließen des Tauschs von Nadelsitz und Nadeleinheit (▶ Seite 258) fort.

7.9.4.3 Abschließen des Tauschs von Nadelsitz und Nadeleinheit

1. Bauen Sie die Halterung der Isolierabdeckung wieder ein, indem Sie die Halterung in die Auflageschienen drücken bis sie einrastet.



Abbildung 157: Einbauen der Halterung der Isolierabdeckung

- 2. Ziehen Sie die Nadeleinheit mit der Hand leicht nach vorn heraus.
- 3. Sichern Sie die Nadeleinheit.
- Wenn die Probenschleife entfernt wurde: Bauen Sie die Probenschleife wieder ein. Folgen Sie den Installationsschritten in Abschnitt Einbauen der Probenschleife (▶ Seite 244).
- 5. Führen Sie die Kapillare von der Probenschleife durch die Halterung der Isolierabdeckung zum Injektionsventil.
- 6. Entsichern Sie die Nadeleinheit.
- 7. Bauen Sie die Isolierabdeckung wieder ein.
- 8. Schalten Sie den Autosampler über den Hauptnetzschalter wieder ein. Der Autosampler führt einen Selbsttest durch. Die Nadel wird initialisiert und justiert ihre Position.
- Abhängig vom Tausch: Stellen Sie die entsprechenden Parameter im Chromatographie-Datensystem ein, zum Beispiel NeedleChanged und NeedleSeatChanged.

7.10 Wartung des Kopfes der Dosiereinheit

7.10.1 Komponenten des Kopfes der Dosiereinheit



Abbildung 158: Übersicht über den Kopf der Dosiereinheit

Nr.	Beschreibung
1	Kopf der Dosiereinheit
2	Befestigungsschrauben
3	Hinterspülauslass
4	Kapillare zum Drucksensor
5	Kapillare zum Injektionsventil
6	Hinterspüleinlass

7.10.2 Entfernen und Anschließen von Kapillaren am Kopf der Dosiereinheit

- 1. Bereiten Sie den Autosampler für die Wartung vor:
 - Schalten Sie den Pumpenfluss ab. Warten Sie, bis das System druckfrei ist, bevor Sie mit der Außerbetriebnahme des Autosamplers fortfahren.
 - Um die Drainagepumpe zu entleeren, lassen Sie die Drainagepumpe laufen, bis im Ablaufschlauch der Drainagepumpe kein Kondenswasser mehr vorhanden ist. Schalten Sie danach die Drainagepumpe aus und entfernen Sie den Peristaltikschlauch aus der Drainagepumpe. Flüssigkeit, die sich im Schlauch der Drainagepumpe befindet, läuft in den Abfall ab.
 - ◆ Fahren Sie die Nadel in die Service-Position und sichern Sie die Nadeleinheit. Siehe Sichern der Nadeleinheit (▶ Seite 248).

Entfernen von Kapillaren vom Kopf der Dosiereinheit

- 1. Entfernen Sie die Injektionsventil-Kapillare vom vorderen Port des Kopfes der Dosiereinheit.
- 2. Entfernen Sie die Drucksensor-Kapillare vom unteren Port des Kopfes der Dosiereinheit.
- 3. Wenn die Kapillaren wieder angeschlossen werden sollen, schützen Sie offene Kapillar-Fittinge mit Schutzkappen.

Anschließen der Kapillaren am Kopf der Dosiereinheit

- 1. Schließen Sie die Injektionsventil-Kapillare an den vorderen Port des Kopfes der Dosiereinheit an.
- Schließen Sie die Drucksensor-Kapillare an den unteren Port des Kopfes der Dosiereinheit an.
 Die Befestigungsschraube des Drucksensors kann für den leichteren Zugang gelöst und wieder festgeschraubt werden.
- 3. Schließen Sie die Wartung des Autosamplers ab:
 - Entsichern Sie die Nadeleinheit. Siehe Entsichern der Nadeleinheit (▶ Seite 249).
 - Starten Sie den Autosampler neu.
 - Bereiten Sie den Autosampler f
 ür den Betrieb vor. Siehe Vorbereiten des Systems f
 ür den Betrieb (▶ Seite 160).

7.10.3 Tauschen des Kopfes der Dosiereinheit

Wann

Kopf der Dosiereinheit ist beschädigt

Erforderliche Teile

Ersatz-Kopf der Dosiereinheit

Erforderliche Werkzeuge

- Schraubendreher, Torx T30
- Pinzette (optional)

- 1. Um den Kopf der Dosiereinheit abzukoppeln, führen Sie das Skript aus: **C22 Kopf der Dosiereinheit auswechseln**.
- 2. Bereiten Sie den Autosampler für die Wartung vor:
 - Schalten Sie den Pumpenfluss ab. Warten Sie, bis das System druckfrei ist, bevor Sie mit der Außerbetriebnahme des Autosamplers fortfahren.
 - Um die Drainagepumpe zu entleeren, lassen Sie die Drainagepumpe laufen, bis im Ablaufschlauch der Drainagepumpe kein Kondenswasser mehr vorhanden ist. Schalten Sie danach die Drainagepumpe aus und entfernen Sie den Peristaltikschlauch aus der Drainagepumpe. Flüssigkeit, die sich im Schlauch der Drainagepumpe befindet, läuft in den Abfall ab.
 - Fahren Sie die Nadel in die Service-Position und sichern Sie die Nadeleinheit. Siehe Sichern der Nadeleinheit (> Seite 248).
- 3. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.
- Entfernen Sie die Kapillaren vom Kopf der Dosiereinheit. Siehe Entfernen und Anschließen von Kapillaren am Kopf der Dosiereinheit (▶ Seite 260).
- Stecken Sie die Hinterspülschläuche vom Kopf der Dosiereinheit ab. Um das Entfernen zu erleichtern, ziehen Sie die Schlauchverbinder mit den Schläuchen vom Kopf der Dosiereinheit unter Verwendung einer Pinzette ab.

Entfernen des Kopfes der Dosiereinheit

- 1. Lösen Sie mit dem Schraubendreher die zwei Schrauben, die den Kopf der Dosiereinheit am Autosampler befestigen.
- 2. Entfernen Sie den Kopf der Dosiereinheit:
 - a) Greifen Sie den Kopf der Dosiereinheit an der Vorderseite. Bewegen Sie den Kopf der Dosiereinheit vorsichtig einige Male nach links und rechts, um den Kolben von der Dosiereinheit zu lösen.
 - b) Überprüfen Sie, ob der Kopf der Dosiereinheit circa 5 mm von der Frontplatte des Autosamplers heraussteht.
 - c) Drücken Sie die Vorderseite des Kopfes der Dosiereinheit vorsichtig nach unten, und ziehen Sie den Kopf von der Frontplatte weg.
 - d) Achten Sie darauf, dass der Kopf der Dosiereinheit mit dem Kolben entfernt wurde.



Abbildung 159: Lösen des Kopfes der Dosiereinheit

Nr.	Beschreibung
1	Kolben
2	Kopf der Dosiereinheit

TIPP Wenn der Kolben in der Dosiereinheit feststeckt, entfernen Sie zuerst den Kolben, bevor Sie mit der Tauschprozedur fortfahren. Siehe Abschnitt .

Einbauen des Kopfes der Dosiereinheit

- 1. Entpacken Sie den Ersatz-Kopf der Dosiereinheit. Lassen Sie die Verschlussfittinge angeschlossen und die Hinterspülschläuche am Kopf der Dosiereinheit miteinander verbunden.
- 2. Setzen Sie den Kopf der Dosiereinheit in die Öffnung an der Innenwand.
- Ziehen Sie die zwei Schrauben mit dem Schraubendreher (Torx T30) am Kopf der Dosiereinheit handfest an bis sie nicht weiter angezogen werden können.
- 4. Entfernen Sie die Verschlussstopfen vom Kopf der Dosiereinheit.
- 5. Stecken Sie die miteinander verbundenen Hinterspülschläuche vom Schlauchverbinder in der Mitte ab. Die Hinterspülschläuche können am Kopf der Dosiereinheit angeschlossen bleiben.
- Schließen Sie die Hinterspülschläuche des neuen Kopfes der Dosiereinheit an die Hinterspülung an. Folgen Sie den Installationsschritten entsprechend für die Hinterspülschläuche, die angeschlossen werden sollen, in Tauschen der Hinterspülschläuche.
- Schließen Sie die Kapillaren an den Kopf der Dosiereinheit an. Siehe Entfernen und Anschließen von Kapillaren am Kopf der Dosiereinheit (> Seite 260).
- 8. Schließen Sie die Wartung des Autosamplers ab:
 - Entsichern Sie die Nadeleinheit. Siehe Entsichern der Nadeleinheit (▶ Seite 249).
 - Starten Sie den Autosampler neu.
 - Bereiten Sie den Autosampler f
 ür den Betrieb vor. Siehe Vorbereiten des Systems f
 ür den Betrieb (
 Seite 160).
- Führen Sie das Skript aus: C22 Kopf der Dosiereinheit auswechseln. Der Autosampler bereitet den Kopf der Dosiereinheit für den Betrieb vor.
- 10. Testen Sie die Hinterspülung auf Dichtigkeit. Siehe Testen der Hinterspülung auf Undichtigkeit (▶ Seite 279).
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

7.11 Wartung der Drainagepumpe

Die Drainagepumpe (Peristaltikpumpe) im Autosampler entfernt jegliche Kondensationsflüssigkeit aus dem Probenraum, die gegebenenfalls während der Probenkühlung entsteht und führt sie durch einen separaten Drainage-Ablauf in den Abfall.



Abbildung 160: Drainagepumpe

Nr.	Beschreibung
1	Hebel der Pumpe
2	Rotor der Drainagepumpe
3	Peristaltikschlauch
4	Schlauchhalterung
5	Kondensationsschlauch zum Probenraum
6	Ablaufschlauch zum Drainage-Ablauf

7.11.1 Testen des Drainagepumpen-Schlauchs

Testen Sie den Drainagepumpen-Schlauch auf Blockierungen und Undichtigkeit.

Wann

In regelmäßigen Abständen

Erforderliche Teile

Spritze (12 mL)

- 1. Schalten Sie die Drainagepumpe aus.
- 2. Ziehen Sie die Spritze mit Luft auf.

3. Ziehen Sie den Ablaufschlauch und den L-Schlauchverbinder vom Peristaltikschlauch ab.



Abbildung 161: Abziehen des Ablaufschlauchs und des L-Schlauchverbinders

- 4. Führen Sie die Spritze in das offene Ende des Peristaltikschlauchs ein.
- 5. Damit die Luft die Drainage-Verschlauchung leicht passieren kann, drücken Sie den Hebel der Drainagepumpe nach oben und halten Sie ihn mit einer Hand in dieser Position.



Abbildung 162: Drücken des Pumpenhebels nach oben

Testen auf Blockierungen

- 1. Drücken Sie den Spritzenstößel, um Luft in den Schlauch zu spritzen.
- 2. Achten Sie darauf, ob ein Widerstand am Stößel vorhanden ist, wenn Sie den Spritzenstößel drücken:

Beobachtung	Schritte
Widerstand ist vorhanden	Der Peristaltikschlauch ist möglicherweise blockiert. Entfernen Sie die Spritze und tauschen Sie den Peristaltikschlauch aus (siehe Tauschen des Drainagepumpen-Schlauchs () Seite 267)).
Es ist kein Widerstand vorhanden	Testen Sie den Schlauch auf Dichtigkeit.

Testen auf Dichtigkeit

1. Ziehen Sie die Spritze mit Luft auf.

2. Ziehen Sie den Kondensationsschlauch und den L-Schlauchverbinder vom Peristaltikschlauch ab.



Abbildung 163: Abziehen des Kondensationsschlauchs und des L-Schlauchverbinders

- 3. Halten Sie ein Ende des Peristaltikschlauchs mit einem Finger zu.
- 4. Drücken Sie den Spritzenstößel, um Luft in den Schlauch zu spritzen.
- 5. Achten Sie darauf, ob sich Druck aufbaut, wenn Sie den Spritzenstößel drücken:

Beobachtung	Schritte
Es ist kein Druck vorhanden	Der Peristaltikschlauch ist möglicherweise undicht oder beschädigt. Entfernen Sie die Spritze und tauschen Sie den Peristaltikschlauch aus (siehe Tauschen des Drainagepumpen- Schlauchs () Seite 267)).
Druck ist vorhanden	Entfernen Sie die Spritze und schließen Sie die L-Schlauchverbinder, den Kondensationsschlauch und den Ablaufschlauch wieder an. Bewegen Sie den Hebel in die Ausgangsstellung zurück.

7.11.2 Tauschen des Drainagepumpen-Schlauchs

Wann

- Alle sechs Monate als präventive Wartungsmaßnahme für die Drainagepumpe
- Der Drainagepumpen-Schlauch ist blockiert (siehe Testen des Drainagepumpen-Schlauchs (▶ Seite 265)).
- Der Drainagepumpen-Schlauch ist undicht oder beschädigt (siehe Testen des Drainagepumpen-Schlauchs (▶ Seite 265)).

Erforderliche Teile

- Ersatz-Peristaltikschlauch für die Drainagepumpe
- 2 L-Schlauchverbinder für den Kondensationsschlauch und den Ablaufschlauch

Erforderliche Werkzeuge

Kapillarschneider (optional)

Vorbereitungen

- 1. Schalten Sie die Drainagepumpe aus.
- Ziehen Sie den Ablaufschlauch und den Kondensationsschlauch von den Schlauchverbindern ab, die sich am Drainagepumpen-Schlauch (Peristaltikschlauch) befinden. Lassen Sie die Schlauchverbinder an den Peristaltikschlauch angesteckt.



Abbildung 164: Abziehen des Ablaufschlauchs und des Kondensationsschlauchs

- 3. Entpacken Sie den Ersatz-Peristaltikschlauch und die L-Schlauchverbinder.
- 4. Schließen Sie die L-Schlauchverbinder an beiden Enden des Ersatz-Peristaltikschlauchs an.

Gehen Sie wie folgt vor

1. Drücken Sie den Hebel der Drainagepumpe nach oben und halten Sie ihn mit einer Hand in dieser Position.



Abbildung 165: Drücken des Pumpenhebels nach oben

- Wickeln Sie mit der anderen Hand den eingebauten Peristaltikschlauch vom Rotor der Drainagepumpe und entfernen Sie ihn von der Drainagepumpe. Bewegen Sie den Hebel in die Ausgangsstellung zurück.
- Prüfen Sie, dass der Ersatz-Peristaltikschlauch die gleiche Länge wie der vorherige Schlauch hat. Schneiden Sie den Ersatz-Peristaltikschlauch falls erforderlich mit dem Kapillarschneider auf die Länge zu.
- Drücken Sie den Ersatz-Peristaltikschlauch in die Schlauchhalterungen. Prüfen Sie, dass der Schlauch korrekt in die Schlauchhalterungen eingelegt ist.
- 5. Drücken Sie den Hebel der Drainagepumpe nach oben. Legen Sie den Ersatz-Peristaltikschlauch zwischen Hebel und Rotor der Drainagepumpe ein und um den Rotor herum.
- 6. Bewegen Sie den Hebel in die Ausgangsstellung zurück.
- 7. Schließen Sie den Ablaufschlauch und den Kondensationsschlauch an die Schlauchverbinder am Peristaltikschlauch an.



Abbildung 166: Anschließen des Ablaufschlauchs und des Kondensationsschlauchs

8. Schalten Sie die Drainagepumpe ein, um sie auf korrekte Funktionsfähigkeit zu prüfen.

7.12 Wartung des Nadel-Waschsystems

Eine Übersicht der Waschschläuche finden Sie unter Nadel-Waschsystem (Übersicht) (▶ Seite 58).

7.12.1 Tauschen der Waschschläuche vom Behälter zum Eingang der Nadel-Waschpumpe

Tauschen Sie die beiden Schläuche zwischen den Behältern für Nadel-Waschflüssigkeit und den Einlass-Ports der Nadel-Waschpumpe im Autosampler wie nachfolgend beschrieben.

Wann

Waschschlauch ist verstopft oder beschädigt

Erforderliche Teile und zusätzliche Utensilien

- Frische Nadel-Waschflüssigkeit
- Ersatz-Waschschläuche mit Beschriftung SWP (rot) und WWP (grün)

Erforderliche Werkzeuge

Kapillarschneider (optional)

TIPP Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.

- 1. Bereiten Sie eine frische Nadel-Waschflüssigkeit vor.
- 2. Entleeren Sie das Waschsystem für das Waschen des Nadeläußeren:
 - a) Schrauben Sie die Deckel der Behälter für die Nadelwaschflüssigkeit ab.
 - b) Entfernen Sie die Waschschläuche (Beschriftung WWP und SWP) zusammen mit dem Deckel aus dem Behälter.
 - Wenn Sie die Schläuche aus den Schlauchführungen entfernen, achten Sie darauf, nicht an den anderen Schläuchen in den Führungen zu ziehen.
 - d) Führen Sie die Waschschläuche in einen Abfallbehälter, um die Schläuche zu entleeren.

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Schlauchführung von den Waschschläuchen.
- 2. Ziehen Sie die Schläuche aus den Deckeln der Behälter.
- 3. Stecken Sie die Waschschläuche von den Einlass-Ports an der Nadel-Waschpumpe ab.



Abbildung 167: Entfernen der Waschschläuche vom Einlass-Port der Nadel-Waschpumpe

- 4. Schließen Sie die Ersatz-Waschschläuche an die Einlass-Ports der Nadel-Waschpumpe an:
 - a) Drücken Sie den steifen Schlauch in den Verbinder.
 - b) Schieben Sie den Silikonschlauch über den Verbinder.
- Führen Sie die Waschschläuche vom Autosampler durch die Schlauchführungen im Autosampler und allen Modulen, die sich über dem Autosampler im Systemturm befinden, zum Solvent Rack.
- 6. Führen Sie die Waschschläuche durch das Führungsloch im Solvent Rack und drücken Sie sie in eine oder mehrere Schlauchführungen.



Abbildung 168: Führungsloch und Schlauchführungen im Solvent Rack

Nr.	Beschreibung
А	Führungsloch (Solvent Rack)
1-4	Schlauchführungen

- 7. Führen Sie jeden Waschschlauch durch eine Schlauchführung.
- 8. Führen Sie jeden Waschschlauch durch eine Öffnung im Deckel eines der Behälter. Die Schlauchführung verhindert, dass der Lösungsmittelschlauch im Behälter verrutscht.
- 9. Verschließen Sie die Öffnungen in den Behälterdeckeln mit den Verschlusskappen.
- 10. Schieben Sie einen Filterhalter mit Adapter auf jeden Waschschlauch.
- Befüllen Sie die Behälter für die Nadel-Waschflüssigkeit mit frischer Nadel-Waschflüssigkeit und setzen Sie das Nadel-Waschsystem wieder in Betrieb. Einzelheiten finden Sie unter Füllen der Behälter für die Nadel-Waschflüssigkeit (> Seite 130).

7.12.2 Tauschen des Waschschlauchs vom Auslass der Nadel-Waschpumpe zum Waschport

Tauschen Sie den Nadel-Waschschlauch zwischen den Auslass-Ports der Nadel-Waschpumpe und dem Waschport wie nachfolgend beschrieben.

Wann

Waschschlauch ist verstopft oder beschädigt

Erforderliche Teile und zusätzliche Utensilien

- Frische Nadel-Waschflüssigkeit
- Ersatz-Waschschlauch

- 1. Bereiten Sie eine frische Nadel-Waschflüssigkeit vor.
- 2. Entleeren Sie das Waschsystem für das Waschen des Nadeläußeren:
 - a) Schrauben Sie die Deckel der Behälter für die Nadelwaschflüssigkeit ab.
 - b) Entfernen Sie die Waschschläuche (Beschriftung WWP und SWP) zusammen mit dem Deckel aus dem Behälter.
 - c) Wenn Sie die Schläuche aus den Schlauchführungen entfernen, achten Sie darauf, nicht an den anderen Schläuchen in den Führungen zu ziehen.
 - d) Führen Sie die Waschschläuche in einen Abfallbehälter, um die Schläuche zu entleeren.

Gehen Sie wie folgt vor

1. Stecken Sie die beiden Enden des Waschschlauchs von den Auslass-Ports der Nadel-Waschpumpe ab.



Abbildung 169: Entfernen der Waschschläuche vom Auslass-Port der Nadel-Waschpumpe

2. Entfernen Sie den Waschschlauch vom Einlass des Waschports.



Abbildung 170: Entfernen des Waschschlauchs vom Waschport

- 3. Schließen Sie den Ersatz-Waschschlauch an die Auslass-Ports der Nadel-Waschpumpe und den Einlass des Waschports an.
- Befüllen Sie die Behälter für die Nadel-Waschflüssigkeit mit frischer Nadel-Waschflüssigkeit und setzen Sie das Nadel-Waschsystem wieder in Betrieb. Einzelheiten finden Sie unter Füllen der Behälter für die Nadel-Waschflüssigkeit (▶ Seite 130).

7.12.3 Tauschen der Waschschläuche vom Behälter zu den Einlass-Ports am Degaser

Tauschen Sie die beiden Schläuche zwischen den Behältern für Nadel-Waschflüssigkeit und den Einlass-Ports des Degasers in der Pumpe wie nachfolgend beschrieben.

Wann

Waschschlauch vom Behälter zum Degaser ist verstopft oder beschädigt

Erforderliche Teile und zusätzliche Utensilien

- Frische Nadel-Waschflüssigkeit
- Ersatz-Waschschläuche mit Beschriftung W (grün) und S (rot)

Erforderliche Werkzeuge

Kapillarschneider (optional)

TIPP Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.

Vorbereitungen

- 1. Bereiten Sie eine frische Nadel-Waschflüssigkeit vor.
- 2. Entleeren Sie das Nadel-Waschsystem:
 - a) Schrauben Sie die Deckel der Behälter für die Nadelwaschflüssigkeit ab.
 - b) Entfernen Sie die Waschschläuche (Beschriftung W und S) zusammen mit dem Deckel aus den Behältern.
 - c) Führen Sie die Waschschläuche in einen Abfallbehälter, um die Schläuche zu entleeren.

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Schlauchführung von den Waschschläuchen.
- 2. Ziehen Sie die Schläuche aus den Deckeln der Behälter.
- 3. Entfernen Sie die Waschschläuche vom Degaser.



4. Schließen Sie die Ersatz-Waschschläuche an die Einlass-Ports des Degasers an.



Nr.	Beschreibung
1	Einlass für Waschschlauch mit Beschriftung ${f W}$ (grün) zum Behälter
2	Einlass für Waschschlauch mit Beschriftung S (rot) zum Behälter
3	Auslass für Waschschlauch mit Beschriftung ${f S}$ (rot) zum Autosampler
4	Auslass für Waschschlauch mit Beschriftung ${f W}$ (grün) zum Autosampler

- 5. Führen Sie die Waschschläuche vom Autosampler durch die Schlauchführungen im Autosampler und allen Modulen, die sich über dem Autosampler im Systemturm befinden, zum Solvent Rack.
- 6. Führen Sie die Waschschläuche durch das Führungsloch im Solvent Rack und drücken Sie sie in eine oder mehrere Schlauchführungen.



Abbildung 172: Führungsloch und Schlauchführungen im Solvent Rack

Nr.	Beschreibung
А	Führungsloch (Solvent Rack)
1-4	Schlauchführungen

- 7. Führen Sie jeden Waschschlauch durch eine Schlauchführung.
- 8. Führen Sie jeden Waschschlauch durch eine Öffnung im Deckel eines der Behälter. Die Schlauchführung verhindert, dass der Lösungsmittelschlauch im Behälter verrutscht.
- 9. Verschließen Sie die Öffnungen in den Behälterdeckeln mit den Verschlusskappen.
- 10. Schieben Sie einen Filterhalter mit Adapter auf jeden Waschschlauch.
- Befüllen Sie die Behälter für die Nadel-Waschflüssigkeit mit frischer Nadel-Waschflüssigkeit und setzen Sie das Nadel-Waschsystem wieder in Betrieb. Siehe Füllen der Behälter für die Nadel-Waschflüssigkeit () Seite 130).

7.12.4 Tauschen der Waschschläuche vom Degaser-Ausgang zum Umschaltventil

Tauschen Sie die beiden Schläuche zwischen den Auslass-Ports des Degasers in der Pumpe und den Einlass-Ports am Umschaltventil des Autosamplers wie nachfolgend beschrieben.

Wann

Waschschlauch ist verstopft oder beschädigt

Erforderliche Teile und zusätzliche Utensilien

- Frische Nadel-Waschflüssigkeit
- Ersatz-Waschschläuche

- 1. Bereiten Sie eine frische Nadel-Waschflüssigkeit vor.
- 2. Entleeren Sie das Nadel-Waschsystem:
 - a) Schrauben Sie die Deckel der Behälter für die Nadelwaschflüssigkeit ab.
 - b) Entfernen Sie die Waschschläuche (Beschriftung W und S) zusammen mit dem Deckel aus den Behältern.
 - c) Führen Sie die Waschschläuche in einen Abfallbehälter, um die Schläuche zu entleeren.

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Schlauchführung von den Waschschläuchen.
- 2. Entfernen Sie die Schläuche vom Umschaltventil des Autosamplers.
- 3. Entfernen Sie die Schläuche vom Auslass-Port des Degasers.
- 4. Schließen Sie die Ersatz-Waschschläuche an die Auslass-Ports des Degasers an.
- 5. Schließen Sie die Waschschläuche an das Umschaltventil des Autosamplers an.
- Befüllen Sie die Behälter für die Nadel-Waschflüssigkeit mit frischer Nadel-Waschflüssigkeit und setzen Sie das Nadel-Waschsystem wieder in Betrieb. Siehe Füllen der Behälter für die Nadel-Waschflüssigkeit () Seite 130).

Sehen Sie dazu auch

■ Nadel-Waschsystem (Übersicht) (▶ Seite 58)

7.12.5 Tauschen des Kugelventils im Nadel-Waschsystem





Nr.	Beschreibung
1	Viper-Kapillare zum Lösungsmittelventil
2	Kugelventil für Nadel-Waschfunktion
3	Umschaltventil für Nadel-Waschfunktion

Wann

Kugelventil ist defekt

Erforderliche Teile

Ersatz-Kugelventil

Gehen Sie wie folgt vor

- 1. Lösen Sie das Viper-Fitting am Kugelventil und entfernen Sie die Viper-Kapillare vom Lösungsmittelventil.
- 2. Schrauben Sie das Kugelventil vom Umschaltventil ab.
- 3. Installieren Sie das Ersatz-Kugelventil.
- 4. Schließen Sie die Viper-Kapillare zum Lösungsmittelventil an. Ziehen Sie die Verbindung fest.
- 5. Purgen Sie den Sampler.

Sehen Sie dazu auch

- Anschlussbelegung der Autosampler-Ventile (> Seite 48)
- Purgen (Übersicht) (> Seite 65)

7.13 Wartung der Hinterspülung

Dieser Abschnitt beschreibt wie Teile der Hinterspülung gewartet und getauscht werden. Folgen Sie den Schritten für die Wartungsprozedur, die Sie ausführen möchten.

7.13.1 Testen der Hinterspülung auf Undichtigkeit

Wann

- In regelmäßigen Abständen
- Wenn der Pumpenkopf bei der Wartung installiert wurde
- Nach Tausch des Kopfes der Dosiereinheit

Erforderliche Utensilien

- Spritze (12 mL)
- Hinterspülflüssigkeit

Gehen Sie wie folgt vor

1. Ziehen Sie den Schlauch vom Hinterspüldetektor ab.



Abbildung 174: Hinterspüldetektor

Nr.	Beschreibung
1	Hinterspüldetektor
2	Schlauchanschluss (Port mit der Kennzeichnung in)

- 2. Führen Sie eine Spritze in das offene Schlauchende ein.
- 3. Damit die Flüssigkeit die Hinterspülung leicht passieren kann, ziehen Sie den Hebel der Hinterspülpumpe nach oben und halten Sie ihn mit einer Hand in dieser Position.
- 4. Ziehen Sie den Spritzenstößel heraus, um Flüssigkeit in die Spritze aufzuziehen.

- 5. Wenn sich circa 10 mL Flüssigkeit in der Spritze befinden, bewegen Sie den Hebel der Hinterspülpumpe in die Ausgangsstellung zurück.
- 6. Drücken Sie den Hinterspülschlauch zum Behälter oberhalb des Kopfs der Dosiereinheit so zusammen, dass dieser Schlauch blockiert wird.
- 7. Drücken Sie den Spritzenstößel leicht, so dass etwas Druck in der Hinterspülung aufgebaut wird.



Abbildung 175: Position, wo der Hinterspülschlauch gehalten werden soll

 Prüfen Sie, ob Flüssigkeit aus den Verbindungen der Hinterspülung, unter den Pumpenköpfen oder unterhalb des Kopfs der Dosiereinheit austritt.

Beobachtung	Abhilfe
Es liegt eine Undichtigkeit an einer Verbindung vor	 Ziehen Sie undichte Verbindungen fest oder tauschen Sie diese aus. Wiederholen Sie den Test.
Es liegt eine Undichtigkeit unter den Pumpenköpfen vor	 Prüfen Sie, dass die Pumpenkopfdichtungen und Hinterspüldichtung korrekt eingebaut sind. Wiederholen Sie den Test.
Es liegt eine Undichtigkeit unterhalb des Kopfs der Dosiereinheit vor	 Prüfen Sie, dass der Kopf der Dosiereinheit korrekt installiert ist. Wenn die Undichtigkeit weiter auftritt, tauschen Sie den Kopf der Dosiereinheit. Wiederholen Sie den Test.
Es liegt keine Undichtigkeit vor	Entfernen Sie die Spritze und schließen den Schlauch wieder am Hinterspüldetektor an (am Port mit der Kennzeichnung in).

Sehen Sie dazu auch

Tauschen des Kopfes der Dosiereinheit (Seite 261)

7.13.2 Tauschen des Hinterspülschlauchs zum Behälter

Tauschen Sie den Spülschlauch zwischen Hinterspülbehälter und Einlass-Port im Autosampler wie nachfolgend beschrieben.



Abbildung 176: Position des Hinterspülschlauchs zum Behälter

Wann

Hinterspülschlauch zum Behälter ist verstopft oder beschädigt

Erforderliche Teile

Silikonschlauch und Schlauchverbinder

Erforderliche Werkzeuge

Kapillarschneider

- 1. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.

Entfernen des Hinterspülschlauchs zum Behälter

1. Ziehen Sie den Hinterspülschlauch mit Schlauchverbinder aus dem Hinterspül-Einlass-Port.



Abbildung 177: Herausziehen des Schlauchverbinders aus dem Hinterspül-Einlass-Port

2. Stecken Sie den Hinterspülschlauch zum Behälter, Schlauchverbinder und Hinterspül-Einlassschlauch voneinander ab. Am Kopf der Dosiereinheit kann der Hinterspülschlauch angeschlossen bleiben.



Abbildung 178: Abstecken des Hinterspülschlauchs vom Hinterspül-Einlassschlauch

- 3. Entfernen Sie den Schlauch vom Hinterspülbehälter:
 - a) Entfernen Sie die Schlauchführung vom Spülschlauch.
 - b) Ziehen Sie den Spülschlauch aus dem Deckel des Behälters.





Nr.	Beschreibung
1	Spülschlauch
2	Schlauchführung
3	Behälterdeckel
4	Verschlusskappen

Verbinden des neuen Hinterspülschlauchs

- 1. Schneiden Sie den Schlauch auf die entsprechende Länge zu, falls notwendig.
- 2. Schließen Sie den Hinterspülschlauch mit einem neuen Schlauchverbinder an den Hinterspül-Einlassschlauch an.
- 3. Drücken Sie den Schlauchverbinder mit den angeschlossenen Schläuchen in den Hinterspül-Einlass-Port.
- Führen Sie den Hinterspülschlauch vom Autosampler durch die Schlauchführungen im Autosampler und allen Modulen, die sich über dem Autosampler im Systemturm befinden, zum Solvent Rack.
- 5. Führen Sie den Hinterspülschlauch durch das Führungsloch im Solvent Rack und befestigen Sie ihn in einer der Schlauchführungen.





Nr.	Beschreibung
А	Führungsloch (Solvent Rack)
1-4	Schlauchführungen

- 6. Schließen Sie den Behälter für die Hinterspülflüssigkeit an den Schlauch an:
 - a) Spülen Sie den Hinterspülbehälter gründlich mit einem hochreinen Lösungsmittel aus.
 - b) Befüllen Sie den Hinterspülbehälter mit frischer Hinterspülflüssigkeit.
 - c) Führen Sie den Hinterspülschlauch durch die Schlauchführung und durch eine Öffnung im Deckel des Behälters für die Hinterspülflüssigkeit. Die Schlauchführung verhindert, dass der Schlauch im Behälter verrutscht. Verschließen Sie die Öffnungen im Behälterdeckel mit Verschlusskappen.

Inbetriebnehmen der Hinterspülung

- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 2. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

Sehen Sie dazu auch

B Hinterspülung (Übersicht) (▶ Seite 60)

7.13.3 Tauschen des Hinterspül-Einlassschlauchs

Tauschen Sie den Spülschlauch zwischen Hinterspül-Einlass-Port im Autosampler und dem Hinterspül-Einlass am Kopf der Dosiereinheit wie nachfolgend beschrieben.



Abbildung 181: Position des Hinterspül-Einlassschlauchs

Wann

Hinterspül-Einlassschlauch ist verstopft oder beschädigt

Erforderliche Teile

Silikonschlauch und Schlauchverbinder

Erforderliche Werkzeuge

Kapillarschneider

Vorbereitungen

- 1. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.

Gehen Sie wie folgt vor

1. Ziehen Sie den Hinterspülschlauch mit Schlauchverbinder aus dem Hinterspül-Einlass oben am Kopf der Dosiereinheit.



Abbildung 182: Entfernen des Hinterspül-Einlassschlauchs am Kopf der Dosiereinheit

2. Ziehen Sie den Hinterspülschlauch mit Schlauchverbinder aus dem Hinterspül-Einlass-Port.



Abbildung 183: Herausziehen des Schlauchverbinders aus dem Hinterspül-Einlass-Port

3. Ziehen Sie den Hinterspülschlauch vom Schlauchverbinder ab. Der Hinterspülschlauch zum Behälter auf der anderen Seite des Schlauchverbinders kann angeschlossen bleiben.



Abbildung 184: Entfernen des Hinterspül-Einlassschlauchs vom Hinterspül-Einlass-Port

- Schließen Sie den neuen Hinterspülschlauch mit einem neuen Schlauchverbinder an den Kopf der Dosiereinheit an. Schneiden Sie den Schlauch auf die entsprechende Länge zu, falls notwendig.
- 5. Führen Sie den Hinterspülschlauch zum Hinterspül-Einlass-Port.
- 6. Schließen Sie den Hinterspül-Einlassschlauch an den Schlauchverbinder an.
- 7. Drücken Sie den Schlauchverbinder mit den angeschlossenen Schläuchen in den Hinterspül-Einlass-Port.

Inbetriebnehmen der Hinterspülung

- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 2. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

7.13.4 Tauschen des Hinterspülpumpen-Einlassschlauchs

Tauschen Sie den Spülschlauch zwischen dem Auslass am Kopf der Dosiereinheit und dem Einlass der Hinterspülpumpe wie nachfolgend beschrieben.



Abbildung 185: Position des Hinterspülpumpen-Einlassschlauchs

Wann

Hinterspülpumpen-Einlassschlauch ist verstopft oder beschädigt

Erforderliche Teile

Silikonschlauch und Schlauchverbinder

Erforderliche Werkzeuge

Kapillarschneider

- 1. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.

Gehen Sie wie folgt vor

1. Ziehen Sie den Schlauchverbinder mit dem Hinterspülpumpen-Einlassschlauch aus dem Hinterspül-Auslass am Kopf der Dosiereinheit.





2. Entfernen Sie den Hinterspülpumpen-Einlassschlauch vom Peristaltikschlauch an der Hinterspülpumpe in der Pumpe.



Abbildung 187: Entfernen des Hinterspülpumpen-Einlassschlauchs

- 3. Schließen Sie einen neuen Schlauchverbinder an den neuen Hinterspülpumpen-Einlassschlauch an. Schneiden Sie den Schlauch auf die entsprechende Länge zu, falls notwendig.
- 4. Schließen Sie den neuen Hinterspülpumpen-Einlassschlauch an den Hinterspülauslass am Kopf der Dosiereinheit an.
- 5. Schließen Sie den Hinterspülpumpen-Einlassschlauch an den Peristaltikschlauch der Hinterspülpumpe an.
Inbetriebnehmen der Hinterspülung

- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 2. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

7.13.5 Tauschen des Peristaltikschlauchs in der Hinterspülpumpe

Tauschen Sie den Peristaltikschlauch in der Hinterspülpumpe wie unten beschrieben.



Abbildung 188: Position des Peristaltikschlauchs in der Hinterspülpumpe

Wann

- Schlauch ist verstopft oder beschädigt
- Alle sechs Monate als präventive Wartungsmaßnahme für die Hinterspülpumpe

Erforderliche Teile

Peristaltikschlauch (PharMed[™]-Schlauch), Schlauchverbinder und -halter

Erforderliche Werkzeuge

Kapillarschneider

Vorbereitungen

- 1. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.

Tauschen des Peristaltikschlauchs



Abbildung 189: Hinterspülpumpe

Nr.	Beschreibung
1	Hebel der Pumpe (angehoben)
2	Rotor der Pumpe
3	Hinterspülschlauch vom Autosampler (Silikonschlauch)
4	Schlauchverbinder
5	Schlauchhalter für Peristaltikschlauch
6	Hinterspülschlauch zum Pumpenkopf (Silikonschlauch)
7	Peristaltikschlauch

- 1. Trennen Sie den Peristaltikschlauch vom Schlauchverbinder, der ihn mit dem Silikonschlauch verbindet.
- 2. Ziehen Sie den Hebel der Hinterspülpumpe nach oben und entfernen Sie den Peristaltikschlauch aus der Pumpe und den Schlauchhalterungen.
- 3. Verbinden Sie den Peristaltikschlauch mit Hilfe der Schlauchverbinder mit dem Silikonschlauch. Schneiden Sie den Schlauch auf die entsprechende Länge zu, falls notwendig.

- 4. Legen Sie den Peristaltikschlauch in die obere rechte Schlauchhalterung ein.
- 5. Ziehen Sie den Hebel der Hinterspülpumpe nach oben und halten Sie ihn mit einer Hand in dieser Position.
- 6. Legen Sie den Schlauch mit der anderen Hand zwischen Hebel und Rotor ein und um den Rotor herum.
- 7. Bewegen Sie den Hebel in die Ausgangsstellung zurück.
- 8. Legen Sie den Peristaltikschlauch in die untere rechte Schlauchhalterung ein.
- 9. Legen Sie den Peristaltikschlauch in die linke Schlauchhalterung ein.
- Prüfen Sie, dass der Peristaltikschlauch korrekt in die Schlauchhalterungen eingelegt ist. Falls nicht, drücken Sie den Schlauch in die Halterung. Vergewissern Sie sich, dass der Schlauch nicht in den Halterungen eingeklemmt oder gequetscht ist.
- 11. Nehmen Sie die Hinterspülung wieder in Betrieb.

Inbetriebnehmen der Hinterspülung

- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 2. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

7.13.6 Tauschen der Hinterspülschläuche zwischen Hinterspülpumpe und Hinterspüldetektor

Tauschen Sie die Hinterspülschläuche zwischen dem Peristaltikschlauch (Hinterspülpumpe) und dem Hinterspüldetektor wie unten beschrieben.





Nr.	Beschreibung
1	Spülschlauch zwischen Peristaltikschlauch und linkem Pumpenkopf
2	Spülschlauch am linken Pumpenkopf
3	Spülschlauch zwischen den Pumpenköpfen
4	Spülschlauch am rechten Pumpenkopf
5	Spülschlauch zwischen rechtem Pumpenkopf und Hinterspüldetektor

Wann

Hinterspülschläuche sind verstopft oder beschädigt

Erforderliche Teile

Silikonschlauch

Erforderliche Werkzeuge

Kapillarschneider

Vorbereitungen

- 1. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.

Gehen Sie wie folgt vor

- 1. Ziehen Sie die Hinterspülschläuche von den Schlauchverbindern ab. Entfernen Sie nicht die Schlauchverbinder.
- 2. Schließen Sie die neuen Hinterspülschläuche an. Schneiden Sie den Schlauch auf die entsprechende Länge zu, falls notwendig.

Inbetriebnehmen der Hinterspülung

- Drehen Sie die Deckel der Lösungsmittelbehälter wieder fest. Die Schlauchführung muss in der Öffnung im Behälterdeckel verbleiben. Ist das nicht der Fall, drücken Sie die Schlauchführung in die Öffnung, so dass der Schlauch im Deckel arretiert wird.
- 2. Prüfen Sie die Schläuche über den gesamten Flussweg hinweg:
 - a) Stellen Sie sicher, dass die Schläuche an keiner Stelle im Flussweg geknickt, eingeklemmt oder gequetscht sind.
 - b) Falls erforderlich, verwenden Sie einen Kapillarschneider, um die Schläuche auf die entsprechende Länge zuzuschneiden. Achten Sie auf einen rechtwinkligen Schnitt zur Längsseite der Leitung.
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

7.13.7 Tauschen des Hinterspüldetektors

Wann

- Verschmutzung der Detektorelektroden
- Eingeschränkte Funktion der Hinterspülung

Erforderliche Teile

Detektoreinheit Hinterspülung, bestehend aus Detektor und Trichter

Vorbereitungen

Drücken Sie zur Vorbereitung der neuen Detektoreinheit Hinterspülung den Hinterspüldetektor in den Trichter. Achten Sie auf die Ausrichtung des Trichters.





Nr.	Beschreibung
1a	Trichter (Unterseite)
1b	Trichter (Oberseite)
2	Hinterspüldetektor

Gehen Sie wie folgt vor





Nr.	Beschreibung
1	Trichter, Hinterspüldetektor
2	Hinterspüldetektor
3	Detektorkabel
4	Hinterspülschlauch

1. Ziehen Sie den Hinterspülschlauch vom Hinterspüldetektor ab (Port mit der Kennzeichnung **in**).

- 2. Ziehen Sie das Detektorkabel vom Port DROP DET ab.
- 3. Fassen Sie die Detektoreinheit Hinterspülung am Trichter und ziehen Sie die Einheit nach oben.
- 4. Drücken Sie die neue Detektoreinheit Hinterspülung in die Pumpe.



Abbildung 193: Einbauen der Detektoreinheit der Hinterspülung

- 5. Schließen Sie den Hinterspülschlauch wieder am Hinterspüldetektor an (Port mit der Kennzeichnung **in**).
- 6. Stecken Sie das Detektorkabel an den Port **DROP DET** an.
- Spülen Sie die Hinterspülung. Führen Sie das Skript aus: B01 -Flüssigkeiten/Lösungsmittel ändern (mit: Hinterspülflüssigkeit hinten und Nur aktualisieren = Aus).

7.14 Prüfen der Autosampler-Ventile auf Dichtigkeit

Wann

- Flüssigkeit tritt an einer Viper-Verbindung aus
- Flüssigkeit tritt an der Ablauföffnung aus



Abbildung 194: Ablauföffnungen an den Autosampler-Ventilen

Erforderliche Werkzeuge

Drehmoment-Schraubendreher mit Steckaufsatz für Viper und Nadelsitz

Gehen Sie wie folgt vor

- Flüssigkeit tritt an der Ablauföffnung eines Ventils aus: Wischen Sie die Flüssigkeit auf und beobachten Sie das Ventil. Treten dann immer noch Flüssigkeitstropfen an der Ablauföffnung auf, wenden Sie sich an den Kundendienst von Thermo Fisher Scientific. Andernfalls fahren Sie mit den nächsten Schritten fort.
- Da Undichtigkeiten in der Regel an Verbindungen auftreten, unterziehen Sie alle Verbindungen im Flussweg einer optischen Prüfung.
- Lösen Sie Kapillarverbindungen am betroffenen Ventil und ziehen Sie sie mit dem Drehmoment-Schraubendreher mit Viper-Steckaufsatz wieder fest.

ACHTUNG

- Lösen und ziehen Sie die Viper-Kapillaren nur mit den Fingern und dem speziellen Drehmoment-Werkzeug fest. Verwenden Sie ausschließlich die für diesen Zweck von Thermo Fisher Scientific zur Verfügung gestellten Werkzeuge. Verwenden Sie das Werkzeug nur für die Verbindung, für die es vorgesehen ist.
- Um Schäden an den Kapillaren und Anschlüssen zu vermeiden, lösen und ziehen Sie die Viper-Kapillaren nur fest, wenn der Systemdruck gleich Null ist.

4. Wenn das Injektionsventil mit dem Nadelsitz das betroffene Ventil ist:

Lösen Sie den Nadelsitz mit dem Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz und ziehen Sie ihn wieder fest.

ACHTUNG—Andere Werkzeuge können den Nadelsitz beschädigen

Der Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz ist zum Festziehen des Nadelsitzes gedacht. Bei Verwendung dieses Werkzeugs wird der Nadelsitz korrekt festgezogen und wird nicht zu fest angezogen. Verwenden Sie kein anderes Werkzeug.

ACHTUNG—Der Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz kann beschädigt werden, wenn er nicht korrekt verwendet wird.

- Setzen Sie den Nadelsitz-Steckaufsatz vollständig in den Drehmoment-Schraubendreher ein.
- Setzen Sie den Drehmoment-Schraubendreher mit Nadelsitz-Steckaufsatz direkt auf den Nadelsitz.
- Drücken oder ziehen Sie das Werkzeug nie mit unnötiger Kraft.
- 5. Ziehen Sie alle Verbindungen im Flussweg mit dem Drehmoment-Schraubendreher mit Viper-Steckaufsatz nach.
- Wenn die Undichtigkeit weiterhin besteht, führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler).
- Flüssigkeitstropfen treten weiterhin am Injektionsventil auf: Tauschen Sie die Nadeleinheit und den Nadelsitz. Siehe Tauschen des Nadelsitzes und der Nadeleinheit (> Seite 251).

7.15 Wartung des Säulenthermostaten

7.15.1 Tauschen der Säulen



VORSICHT—Heiße Säulen

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

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

Wann

Wechseln zwischen Säulen

Erforderliche Teile

Säule, die für Ihre Anwendung erforderlich ist

Vorbereitungen

1. Schalten Sie den Pumpenfluss ab.

Gehen Sie wie folgt vor

- 1. Entfernen Sie die Kapillaren vom Eingang und Ausgang der Säule.
- Entfernen Sie die Säule, die Sie tauschen möchten, aus dem Säulenbefestigungsclip. Drücken Sie den Ring in Richtung der Pfeile, die auf dem Ring angegeben sind, leicht zusammen, um ihn vom Unterteil des Säulenbefestigungsclips abzuziehen.



Abbildung 195: Entfernen des Rings

- 3. Stellen Sie die Kapillarverbindungen an Eingang und Ausgang der neuen Säule her.
- 4. Legen Sie die Säule in das Unterteil des Befestigungsclips und schieben Sie den Ring auf.

5. Führen Sie die Kapillaren durch eine der in der Abbildung gezeigten Kapillarführungen.



Abbildung 196: Kapillarführungen

- Um die S\u00e4uleneigenschaften einzustellen, f\u00fchren Sie die Skripte aus: A03 - Trenns\u00e4ulentyp festlegen und A04 -Trenns\u00e4ulenspezifikationen festlegen (und A05 – Trap-S\u00e4ulenspezifikationen festlegen nach Bedarf).
- 7. Schalten Sie den Pumpenfluss ein.

7.15.2 Visuelle Prüfung der Säulenschaltventile auf Dichtigkeit

Das Vanquish-Ventil und das VICI-Ventil besitzen je eine Ablauföffnung.



Abbildung 197: Ablauföffnungen am Vanquish-Ventil (links) und am VICI-Ventil (rechts)

Wann

- Flüssigkeit tritt an der Ablauföffnung aus
- Flüssigkeit hat sich in der Leckage-Wanne gesammelt

Gehen Sie wie folgt vor

- Da Undichtigkeiten in der Regel an Verbindungen auftreten, unterziehen Sie alle Verbindungen im Flussweg einer optischen Prüfung.
- 2. Entfernen Sie die Kapillaren und schließen Sie die Kapillaren erneut an.
- 3. Ziehen Sie alle Verbindungen im Flussweg nach.
- 4. Prüfen Sie die Ventile visuell auf Undichtigkeit.
- Führen Sie das Skript aus: D02 Systemdichtigkeit testen (was getestet werden soll = Sampler). Dieses Skript testet auch die Säulenschaltventile auf Undichtigkeit.
- Bleibt die Undichtigkeit bestehen, tauschen Sie abhängig vom Ventil das Vanquish-Ventil oder den Statoren und/oder die Rotordichtung des VICI-Ventils.

Für den Tausch des Vanquish-Ventils, siehe Tauschen des Vanquish-Schaltventils (Seite 300).

Für den Tausch des Statoren und/oder der Rotordichtung des VICI-Ventils, siehe Tauschen des Statoren und/oder der Rotordichtung des VICI-Schaltventils (▶ Seite 305).

7.15.3 Tauschen des Vanquish-Schaltventils

Wann

- Schaltventil ist defekt
- Wechseln zwischen verschiedenen Ventiltypen

Erforderliche Teile

- Vanquish-Schaltventil, das für Ihre Anwendung erforderlich ist
- 4 Schrauben

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Vorbereitungen

- 1. Schalten Sie den Pumpenfluss ab.
- 2. Lösen Sie alle Verbindungen am Schaltventil.
- 4. Entfernen Sie die Verschlussfittinge noch nicht von den Ventilports des Ersatz-Schaltventils.

Ausbauen des Vanquish-Schaltventils

Abbildung 198: 4 Schrauben am Schaltventil

2. Ziehen Sie das Schaltventil vorsichtig aus dem Gehäuse. Um das Ventil leichter zu entfernen, schrauben Sie dazu einen Verschlussfitting in einen Ventilport.

1. Lösen Sie die 4 Schrauben am Schaltventil. Bewahren Sie die Schrauben für den Einbau des neuen Schaltventils auf.



Abbildung 199: Ausbauen des Vanquish-Schaltventils



Abbildung 200: Trennen des Kabels vom Stecker

Wenn nach dem Tausch kein Ventil installiert wird

1. Dichten Sie den Kabelstecker mit einem wasserfesten Klebeband ab.



Abbildung 201: Abdichten des Kabelsteckers

- 2. Schieben Sie das Kabel zurück ins Gehäuse.
- 3. Legen Sie die Abdeckung über die Öffnung des Ventilgehäuses.
- 4. Ziehen Sie die 4 Schrauben an der Abdeckung fest.

Installieren des Vanquish-Schaltventils

1. Schließen Sie das neue Schaltventil am Kabel an.

Abbildung 202: Anschließen des Kabels

2. Schieben Sie das Kabel zurück ins Gehäuse.



3. Schieben Sie das Schaltventil vorsichtig ins Gehäuse hinein.

Abbildung 203: Installieren des Vanquish-Schaltventils

- 4. Prüfen Sie die Ausrichtung des Schaltventils: Port 3 muss nach oben zeigen.
- 5. Ziehen Sie die 4 Schrauben am Schaltventil mit einem Schraubendreher fest.
- 6. Entfernen Sie die Verschlussfittinge von den Ventilports.
- 7. Stellen Sie die Verbindungen je nach Erfordernis für Ihre Anwendung her.
- 8. Schalten Sie den Säulenthermostaten ein.
- 9. Prüfen Sie die Ventile visuell auf Undichtigkeit.
- 10. Führen Sie das Skript aus: **D02 Systemdichtigkeit testen (was getestet werden soll = Sampler)**. Dieses Skript testet auch die Säulenschaltventile auf Undichtigkeit.

7.15.4 Tauschen des VICI-Schaltventils

Sie können VICI-Schaltventile nur dort tauschen, wo ein VICI-Ventilantrieb installiert ist. Um den VICI-Ventilantrieb zu installieren, wenden Sie sich an den Kundendienst von Thermo Fisher Scientific. Eine Übersicht des VICI-Schaltventils finden Sie unter Übersicht der Säulenschaltventile (▶ Seite 54).

Voraussetzung

Der VICI-Ventilantrieb und der Adapterrahmen sind installiert (durch Servicetechniker erfolgt).

Wann

- Schaltventil ist defekt
- Wechseln zwischen verschiedenen Ventiltypen

Erforderliche Teile

• VICI-Schaltventil, das für Ihre Anwendung erforderlich ist

Erforderliche Werkzeuge

• Schraubendreher, Torx T10

Vorbereitungen

- 1. Schalten Sie den Pumpenfluss ab.
- 2. Lösen Sie alle Verbindungen am VICI-Schaltventil.
- 3. Entfernen Sie die Verschlussfittinge noch nicht von den Ventilports des Ersatz-Schaltventils.
- 4. Um den installierten Ventilkopf vom Ventilflansch zu entfernen, schrauben Sie einen Verschlussfitting in einen Ventilport und ziehen Sie den Ventilkopf mit Hilfe des Verschlussfittings nach vorn ab.

Wenn nach dem Tausch kein Ventil installiert wird

- 1. Lösen Sie die 4 Schrauben des Adapterrahmens.
- 2. Entfernen Sie den Adapterrahmen.
- 3. Legen Sie die Abdeckung über die Öffnung des Ventilgehäuses.
- 4. Ziehen Sie die 4 Schrauben an der Abdeckung fest.

Installieren des VICI-Schaltventils

1. Richten Sie die beiden Stifte auf der Rückseite des Ventilklemmrings an den entsprechenden Öffnungen im Ventilflansch aus.



Abbildung 204: Ausrichten des Ventilkopfes am Flansch

2. Drücken Sie den Ventilkopf auf den Ventilflansch.

- Um die Achse mit dem Ventilkopf auszurichten, halten Sie je nach Ventilposition die Taste VALVE LEFT oder VALVE RIGHT auf der Tastatur gedrückt. Die Funktion "Catch Valve" wird aktiviert und Sie können die Taste loslassen.
- 4. Wenn die Achse ausgerichtet ist, rastet der Ventilkopf in den Ventilflansch ein.
- 5. Entfernen Sie die Verschlussfittinge von den Ventilports.
- 6. Stellen Sie die Verbindungen je nach Erfordernis für Ihre Anwendung her.
- 7. Konfigurieren Sie das VICI-Schaltventil im Chromatographie-Datensystem.
- 8. Prüfen Sie die Ventile visuell auf Undichtigkeit.
- Führen Sie das Skript aus: D02 Systemdichtigkeit testen (was getestet werden soll = Sampler). Dieses Skript testet auch die Säulenschaltventile auf Undichtigkeit.

7.15.5 Tauschen des Statoren und/oder der Rotordichtung des VICI-Schaltventils

Wann

- Schaltventil ist aufgrund eines beschädigten Statoren und/oder einer beschädigten Rotordichtung undicht
- Rotordichtung ist verstopft

Erforderliche Teile

- Ersatz-Stator und/oder Ersatz-Rotordichtung
- Viper-Fitting oder Verschlussfittinge aus Kunststoff

Erforderliche Werkzeuge

Sechskantschlüssel, Größe 9/64"

Vorbereitungen

- 1. Schalten Sie den Pumpenfluss ab.
- 2. Lösen Sie alle Verbindungen am VICI-Schaltventil.
- 3. Schrauben Sie einen Viper-Fitting oder Verschlussfittinge aus Kunststoff in den Ventilkopf, um den installierten Ventilkopf vom Ventilflansch zu entfernen.

4. Ziehen Sie den Ventilkopf mit Hilfe des Viper-Fittings oder der Verschlussfittinge nach vorn ab.

Gehen Sie wie folgt vor

 Entfernen Sie die 2 Statorschrauben mit einem Sechskantschlüssel. Lösen Sie die Schrauben so lange abwechselnd, bis Sie sie entfernen können.



Abbildung 205: Statorschrauben

2. Entfernen Sie den Stator vom Ventilkörper. Legen Sie den Stator auf seiner Außenseite ab, um die Dichtungsfläche nicht zu beschädigen.



Abbildung 206: VICI-Schaltventil

Nr.	Beschreibung
1	Stator
2	Ventilkörper

3. Entfernen Sie die Rotordichtung vorsichtig mit der Hand.

ACHTUNG

Um eine Beschädigung der Dichtungsflächen des Ventils zu vermeiden, verwenden Sie keine spitzen oder scharfkantigen Werkzeuge. Selbst kleinste Kratzer können die Dichtigkeit des Ventils beeinträchtigen.

4. Prüfen Sie die Dichtungsflächen der Rotordichtung und des Stators auf mögliche Kratzer. Wenn mit bloßem Auge Kratzer auf den Dichtungsflächen erkennbar sind, tauschen Sie die Rotordichtung und/oder den Stator.

- 5. Setzen Sie die neue Rotordichtung ein:
 - Beachten Sie die Ausrichtung der Dichtung. Die Dichtungsfläche muss mit den Flusskanälen nach außen zeigen. Das Muster ist asymmetrisch, um zu verhindern, dass die Rotordichtung falsch eingesetzt wird.
 - Achten Sie darauf, dass keine Verunreinigungen an den Ventilteilen haften. Selbst kleinste Partikel können zu Schäden am Ventil führen und die Dichtigkeit beeinträchtigen.
- Setzen Sie den Stator wieder auf und ziehen Sie die beiden Schrauben mit jeweils einer 1/4-Umdrehung so lange abwechselnd an, bis sie angezogen sind.

TIPP Ziehen Sie die Schrauben nicht zu fest an. Sie dienen lediglich der Befestigung und haben keinen Einfluss auf die Dichtkraft. Die Abdichtung erfolgt automatisch, wenn der Stator auf den Ventilkörper trifft.

- 7. Installieren Sie den Ventilkopf wie in dem entsprechenden Abschnitt unter Tauschen des VICI-Schaltventils (▶ Seite 303) beschrieben.
- 8. Prüfen Sie die Ventile visuell auf Undichtigkeit.
- Führen Sie das Skript aus: D02 Systemdichtigkeit testen (was getestet werden soll = Sampler). Dieses Skript testet auch die Säulenschaltventile auf Undichtigkeit.
- Wenn die Undichtigkeit weiter auftritt, tauschen Sie das Ventil. Siehe Tauschen des VICI-Schaltventils (▶ Seite 303).

7.15.6 Tauschen des passiven Vorheizers

Wann

Passiver Vorheizer defekt

Erforderliche Teile

Passiver Vorheizer

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Gehen Sie wie folgt vor



VORSICHT—Heiße Oberflächen

Der Säulenthermostat und der Vorheizer können heiß werden. Das Berühren von heißen Komponenten kann zu Verbrennungen führen. Lassen Sie den Säulenthermostaten und den Vorheizer abkühlen, bevor Sie Servicearbeiten ausführen.

- 1. Entfernen Sie die Auslasskapillare des passiven Vorheizers vom Säuleneingang.
- 2. Entfernen Sie die Einlasskapillare des passiven Vorheizers vom Autosampler oder Säulenschaltventil.
- Lösen Sie die zwei Schrauben und Unterlegscheiben am passiven Vorheizer mit dem Schraubendreher und entfernen Sie den passiven Vorheizer.
- 4. Befestigen Sie den neuen passiven Vorheizer im Säulenraum. Siehe Installieren des passiven Vorheizers (▶ Seite 92).

7.16 Wartung des System Controllers und des Displays

Dieser Abschnitt beschreibt wie der System Controller und das Display gewartet werden.

7.16.1 Tauschen des System Controllers

- 1. Entfernen Sie den System Controller. Siehe Entfernen des System Controllers (▶ Seite 309).
- 2. Installieren Sie den System Controller. Siehe Installieren des System Controllers (▶ Seite 310).
- 3. Schließen Sie den System Controller an. Siehe Anschließen des System Controllers (▶ Seite 312).
- Wenn Sie den System Controller getauscht haben: Konfigurieren Sie Ihr System im Vanquish User Interface erneut. Informationen zu den Konfigurationsskripten finden Sie unter Vorbereiten des Systems für den ersten Betrieb (▶ Seite 138).

7.16.1.1 Entfernen des System Controllers

Gehen Sie wie folgt vor

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

- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (>> Seite 73).
- 2. Schieben Sie den Systemturm in eine geeignete Position.
- 3. Stecken Sie die Signal- und Netzkabel vom System Controller ab.
- 4. Drücken Sie beide Federklemmen gleichzeitig und ziehen Sie den System Controller kräftig heraus.



Abbildung 207: Drücken der Federklemmen auf der linken und rechten Seite des System Controllers

7.16.1.2 Installieren des System Controllers

Erforderliche Teile

VSC-Ein- und Ausschalter

Erforderliche Werkzeuge

- Zange
- Optional: Schraubendreher, Torx T10

Gehen Sie wie folgt vor

Der Steckplatz für den System Controller befindet sich auf der Rückseite des Systemsockels.

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

- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (>> Seite 73).
- 2. Schieben Sie den Systemturm in eine geeignete Position.
- 3. Stecken Sie die SD-Karte hinein und beachten Sie Folgendes:
 - Korrekte Ausrichtung der Kerbe (siehe Abbildung)
 - Stecken Sie die SD-Karte gerade in das hinter der Aussparung liegende Innengehäuse, bevor Sie die Karte ganz einschieben.



Abbildung 208: Einstecken der SD-Karte

4. Heben Sie den System Controller leicht an und schieben Sie ihn auf der Rückseite des Systemturms mit Kraft in den Systemsockel bis der System Controller mit einem Klick einrastet.



Abbildung 209: Einschieben des System Controllers

5. *Falls der System Controller trotz kräftigem Druck nicht einrastet*: Entfernen Sie die Halteklammer durch Lösen der 3 Schrauben mit dem Schraubendreher.



Abbildung 210: Entfernen der Halteklammer

Nr.	Beschreibung
1	Halteklammer
2	Schrauben

ACHTUNG—Beschädigung des System Controllers

Ohne Halteklammer kann der System Controller beim Anheben des Systems herunterfallen. Dadurch könnte der System Controller beschädigt werden.

- 6. Schließen Sie die Schnittstellenkabel an. Siehe Anschließen des System Controllers (▶ Seite 312).
- Wenn Sie den System Controller getauscht haben: Konfigurieren Sie Ihr System im Vanquish User Interface erneut. Informationen zu den Konfigurationsskripten finden Sie unter Vorbereiten des Systems für den ersten Betrieb (> Seite 138).

7.16.1.3 Anschließen des System Controllers

Folgende Anschlüsse befinden sich auf der kurzen Seite des System Controllers:



Abbildung 211: Anschlüsse auf der kurzen Seite

Nr.	Beschreibung
1	Typenschild
2	Vanquish System Controller (VSC)-Port
3	System Interlink-Port Für den Anschluss an den Systemsockel und an ein Modul im HPLC-System
4	USB-Port (Universal Serial Bus) (Stecker Typ "A") Deaktiviert - Nicht verwenden
5	Local Area Network (LAN)-Port Für den Anschluss an ein firmeninternes Netzwerk.
6	Hauptnetzschalter (Ein/Aus)
7	Netzbuchse

Die folgenden Anschlüsse und Steckplätze befinden sich auf der langen Seite des System Controllers:



Abbildung 212: Anschlüsse auf der langen Seite

Nr.	Beschreibung
1	Netzbuchse für das Display
2	Datensignal-Port für das Display
3	System Interlink-Ports Deaktiviert
4	USB-Port (Stecker Typ "A") Deaktiviert - Nicht verwenden
5	Schild mit Angabe von Teilenummer, Modulnamen und Revisionsnummer.
6	Storage Device (SD)-Kartensteckplatz

ACHTUNG

- Verwenden Sie keine beschädigten Kommunikationskabel. Tauschen Sie das Kabel aus, wenn Sie den Verdacht haben, das Kabel sei beschädigt.
- Verwenden Sie nur die von Thermo Fisher Scientific f
 ür den Anschluss des Systems angebotenen Anschlusskabel, um eine einwandfreie Funktion der Verbindung zu gew
 ährleisten.

Erforderliche Teile

- Netzkabel
- Kurzes VSC-Kabel (Länge: 8 cm)
- System-Interlink-Kabel (Länge: 50 cm)
- Wenn Sie von Ihrem firmeninternen Netzwerk aus auf den System Controller zugreifen wollen: LAN-Kabel (Länge: 5 m)

Gehen Sie wie folgt vor

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 das Gerät über den Netzschalter ausgeschaltet ist.
- 2. Verbinden Sie das Netzkabel mit der Netzbuchse am Gerät.
- 3. Verbinden Sie das andere Ende des Netzkabels mit einer geeigneten Stromquelle.
- 4. Schließen das kurze VSC-Kabel an den **VSC**-Port des Systemsockels und an den **VSC**-Port des System Controllers an.



Abbildung 213: Anschließen der VSC- und System-Interlink-Kabel

ACHTUNG—Schäden am Vanquish-Gerät

Die **System Interlink**- und die **LAN**-Ports besitzen denselben Steckertyp. Wenn Sie zum Beispiel ein LAN-Kabel an einen **System Interlink**-Port eines Vanquish-Moduls anschließen, kann dies das Vanquish-Modul und den System Controller beschädigen.

Verbinden Sie nur Ports desselben Schnittstellentyps.

 Schließen das System-Interlink-Kabel an den System Interlink-Port des Systemsockels und an den System Interlink-Port auf der kurzen Seite des System Controllers an.

- 6. Verbinden Sie den anderen System Interlink-Port auf der kurzen Seite des System Controllers mit dem System Interlink-Port der Pumpe.
- 7. Schließen Sie ein LAN-Kabel an den LAN-Port des System Controllers an.
- Schließen Sie die andere Seite des LAN-Kabels an den LAN-Port der Netzwerkkarte im Datensystemrechner, auf dem die Chromatographie-Software installiert ist.
- 9. Schieben Sie den Systemturm zurück in Betriebsposition.
- 10. Sperren Sie den Systemsockel mit den Systemsockelschlüsseln (siehe Sperren des Systemsockels (▶ Seite 118)).

7.16.2 Tauschen der SD-Karte

TIPP Verwenden Sie nur bei Thermo Fisher Scientific bestellte SD-Karten, um die Kompatibilität der SD-Karte mit dem System Controller sicherzustellen.

Wann

Eine Meldung auf dem Vanquish User Interface fordert Sie auf, die SD-Karte zu tauschen.

Erforderliche Utensilien

- Neue SD-Karte
- Optional: Stift oder Schlitzschraubendreher

Vorbereitungen

- 1. Booten Sie den System Controller und warten Sie, bis die Status-LED des System Controllers leuchtet.
- 2. Prüfen Sie, ob die Meldung erneut erscheint. Wenn ja, fahren Sie mit dem nächsten Schritt fort.
- 3. Schalten Sie den System Controller über den Ein-/Ausschalter des System Controllers aus.

Gehen Sie wie folgt vor

- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (> Seite 73).
- 2. Schieben Sie den Systemturm in eine geeignete Position.

3. Drücken Sie die SD-Karte ein wenig ein, bis die SD-Karte ausgegeben wird. Verwenden Sie hierzu einen Stift oder Schlitzschraubendreher, wenn nötig.



Abbildung 214: Entfernen der SD-Karte

- 4. Entfernen Sie die alte SD-Karte.
- 5. Stecken Sie die SD-Karte hinein und beachten Sie Folgendes:
 - Korrekte Ausrichtung der Kerbe (siehe Abbildung)
 - Stecken Sie die SD-Karte gerade in das hinter der Aussparung liegende Innengehäuse, bevor Sie die Karte ganz einschieben.



Abbildung 215: Einstecken der SD-Karte

- 6. Schieben Sie den Systemturm zurück in Betriebsposition.
- 7. Sperren Sie den Systemsockel mit den Systemsockelschlüsseln (siehe Sperren des Systemsockels (▶ Seite 118)).

7.16.3 Tauschen des Displays

- Bauen Sie das alte Display ab. Siehe Abbauen des Displays (▶ Seite 317).
- Montieren Sie das neue Display. Siehe Montieren des Displays (▶ Seite 100).
- 3. Schließen Sie die Display-Kabel an. Siehe Anschließen der Display-Kabel (▶ Seite 111).

7.16.3.1 Abbauen des Displays

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Gehen Sie wie folgt vor

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

- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (> Seite 73).
- 2. Schieben Sie den Systemturm in eine geeignete Position.
- 3. Ziehen Sie die zwei Display-Stecker vom System Controller ab.
- 4. Entfernen Sie die Kabelabdeckung.
- 5. Entfernen Sie das Display-Kabel aus der Schiene und aus den Kabelhalterungen.
- 6. Halten Sie den Nutenstein mit der rechten Hand und von außen entfernen Sie die Schraube aus einer der Bohrungen, die mit einem Kreis gekennzeichneten sind, mit der linken Hand.



Abbildung 216: Lösen der Nutensteinschraube

- 7. Entfernen Sie den Nutenstein.
- 8. Lösen Sie die 2 Schrauben in den zwei oberen Bohrungen.



Abbildung 217: Lösen der Schrauben

- 9. Schieben Sie das Display zur System-Rückseite und entfernen Sie es.
- 10. Befestigen Sie die oberen 2 Schrauben und befestigen Sie den Nutenstein mit einer Schraube.



Abbildung 218: Die am Display zu befestigenden Schrauben

7.17 Tauschen des Solvent Racks

Erforderliche Teile

- Ersatz-Solvent Rack
- Fusselfreies Tuch

Erforderliche Werkzeuge

Schraubendreher, Torx T10

Vorbereitungen

- 1. Schalten Sie das System über den System-Ein-/Ausschalter (unterer Schalter am Systemsockel) aus.
- 2. Schalten Sie den System-Controller über seinen Ein-/Ausschalter (oberer Schalter am Systemsockel) aus.
- 3. Entleeren Sie die Hinterspülung:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
 - c) Um den Hinterspülschlauch leerlaufen zu lassen, ziehen Sie den Hebel der Hinterspülpumpe nach oben. Bewegen Sie den Hebel in die Ausgangsstellung zurück, wenn der Schlauch leer ist.
- 4. Entfernen und leeren Sie die Lösungsmittelschläuche:
 - a) Schrauben Sie die Deckel der Lösungsmittelbehälter ab.
 - b) Entfernen Sie die Lösungsmittelschläuche zusammen mit den Deckeln aus den Behältern.
 - c) Schützen Sie die Lösungsmittelfilter in geeigneter Form gegen Verschmutzung.
 - d) Entfernen Sie die Lösungsmittelschläuche von den Pumpeneingängen und legen Sie die freien Enden der Lösungsmittelschläuche in die Leckage-Wanne, um die Schläuche zu entleeren.
 - e) Verschließen Sie die Anschlüsse für Lösungsmittel an der Pumpe mit geeigneten Verschlussstopfen.
- 5. Nehmen Sie Flüssigkeit, die aus den Schläuchen austritt, mit einem fusselfreien Tuch auf.
- 6. Entfernen Sie die Flüssigkeitsbehälter aus dem Solvent Rack.

Gehen Sie wie folgt vor

- 1. Lösen Sie mit einem Schraubendreher die Schrauben, die die Blechklammer am Systemmodul unterhalb des Solvent Racks befestigen.
- 2. Entfernen Sie die Blechklammer.

TIPP Bewahren Sie die Blechklammer und die Schrauben auf. Sie benötigen diese gegebenenfalls bei einem künftigen Transport.



Abbildung 219: Entfernen der Blechklammer (hier: Autosampler)

- 3. Heben Sie das Solvent Rack vorne an und ziehen Sie das Solvent Rack nach vorne heraus.
- 4. Ziehen Sie die Schläuche aus dem Schlauchkanal des Solvent Racks heraus.
- 6. Schieben Sie die Schläuche durch den Schlauchkanal des neuen Solvent Racks.
- 7. Setzen Sie das neue Solvent Rack ca. 5 cm vor dem Ende der Schiene auf dem obersten Systemmodul auf. Schieben Sie das Solvent Rack nach hinten, bis es einrastet.
- 8. Setzen Sie den Einsatz für Flaschen in das Solvent Rack ein (siehe Einsetzen des Einsatzes für Flaschen (▶ Seite 323)), wenn vorhanden.
- 9. Stellen Sie die Flüssigkeitsbehälter auf das Solvent Rack.
- 10. Verbinden Sie die Schläuche mit den Flüssigkeitsbehältern.
- 11. Verbinden Sie die Lösungsmittelschläuche mit den Pumpeneingängen.
- 12. Schalten Sie den System-Controller über seinen Ein-/Ausschalter (oberer Schalter am Systemsockel) ein.

- 13. Schalten Sie das System über den System-Ein-/Ausschalter (unterer Schalter am Systemsockel) ein.
- 14. Bereiten Sie das System für die Erstinbetriebnahme vor. Siehe Vorbereiten des Systems für den ersten Betrieb (▶ Seite 138).

7.18 Tauschen des Einsatzes für Flaschen

7.18.1 Entfernen des Einsatzes für Flaschen

Erforderliche Teile

Fusselfreies Tuch

Gehen Sie wie folgt vor

- 1. Schalten Sie das System über den System-Ein-/Ausschalter (unterer Schalter am Systemsockel) aus.
- 2. Schalten Sie den System-Controller über seinen Ein-/Ausschalter (oberer Schalter am Systemsockel) aus.
- 3. Entfernen Sie die Hinterspülschläuche:
 - a) Schrauben Sie den Deckel des Behälters für die Hinterspülflüssigkeit ab.
 - b) Entfernen Sie den Hinterspülschlauch zusammen mit dem Deckel vom Behälter für die Hinterspülflüssigkeit.
- 4. Entfernen Sie die Lösungsmittelschläuche:
 - a) Schrauben Sie die Deckel der Lösungsmittelbehälter ab.
 - b) Entfernen Sie die Lösungsmittelschläuche zusammen mit den Deckeln aus den Behältern.
 - c) Schützen Sie die Lösungsmittelfilter in geeigneter Form gegen Verschmutzung.
- 5. Nehmen Sie Flüssigkeit, die aus den Schläuchen austritt, mit einem fusselfreien Tuch auf.
- 6. Entfernen Sie die Flüssigkeitsbehälter aus dem Solvent Rack.
- 7. Drücken Sie die Flügel des Einsatzes für Flaschen von unten kräftig über die Solvent Rack-Reling und entfernen Sie den Einsatz für Flaschen.



Abbildung 220: Entfernen des Einsatzes für Flaschen

Nr.	Beschreibung
1	Flügel

7.18.2 Einsetzen des Einsatzes für Flaschen

Erforderliche Teile

Einsatz für Flaschen

Gehen Sie wie folgt vor

- 1. Schieben Sie den hinteren und den linken Flügel in den Spalt unter der Reling.
- 2. Drücken Sie den rechten Flügel mit Kraft unter die Reling.



Abbildung 221: Einsetzen des Einsatzes für Flaschen

- 3. Setzen Sie die Flüssigkeitsbehälter in den Einsatz für Flaschen des Solvent-Racks ein.
- 4. Verbinden Sie die Schläuche mit den Flüssigkeitsbehältern.
- 5. Schalten Sie das System und den den System-Controller über die Ein-/Ausschalter vorne links am Systemsockel ein.

7.19 Tauschen der Sicherungen und Türen

7.19.1 Tauschen der Sicherungen

Wann

Durchgebrannte Sicherungen

Erforderliche Teile

Sicherungen (2 Sicherungen, 5 AT, 230 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.





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.

7.19.2 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 Schritten in diesem Abschnitt.

Erforderliche Teile

Ersatztür

Vorbereitungen

Falls sich die zu tauschende Tür direkt unterhalb des Solvent Racks befindet, gehen Sie folgendermaßen vor:

- 1. Entfernen Sie alle Behälter aus dem Solvent Rack.
- 2. Entfernen Sie falls erforderlich die Blechklammer, mit der das Solvent Rack am Autosampler befestigt ist. Siehe Anleitung in Einbau des Säulenthermostaten in den Systemturm (▶ Seite 86).

Falls das Display neben der zu tauschenden Tür installiert ist, lösen Sie gegebenenfalls den Nutenstein des Displays und bewegen Sie das Display nach hinten weg.

Informationen zum Lösen des Nutensteins des Displays finden Sie in Abbauen des Displays (> Seite 317).

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. Falls sich die zu tauschende Tür direkt unterhalb des Solvent Racks befindet, heben Sie die vordere Kante des Solvent Racks leicht an.
- 2. 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 223: Aushängen einer Tür

Nr.	Beschreibung
1	Türangel am Gehäuse
2	Aufnahmenut der Tür

- 3. Kippen Sie die Tür leicht nach außen, weg vom Gehäuse, und entfernen Sie die Tür.
- 4. Zum Einbau der Tür richten Sie die Tür mit den Türangeln am Gehäuse aus. Klemmen Sie keine Schläuche oder Kapillaren zwischen Tür und Gehäuse ein.
- 5. 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.

7.20 Tauschen des Moduleinschubs

Wenn Sie das Gerät an einen anderen Ort transportieren möchten oder versenden müssen, können Sie den Moduleinschub aus dem Gehäuse entfernen. Das Gehäuse verbleibt im Systemturm. Um ein defektes Modul zurückzuschicken, installieren Sie den Moduleinschub in das Gehäuse des Ersatzmoduls.

7.20.1 Vorbereiten für das Entfernen der Pumpe

Allgemeine Vorbereitungen

- 1. Führen Sie das Skript aus: **B05 Außerbetriebnahme (langfristig)** mit dem folgenden Parameter: **Zu entfernendes Modul = Pumpe**.
- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (> Seite 73).
- 3. Schieben Sie den Systemturm in eine geeignete Position.

Außerbetriebnahme

- 1. Schalten Sie *alle* Module über ihren Hauptnetzschalter aus. Das Drücken des System-Ein- /Ausschalters reicht nicht aus, um die Geräte vollständig auszuschalten.
- 2. Ziehen Sie den Netzstecker.
- 3. Entfernen Sie alle Signalkabel zu anderen Geräten.

Flussverbindungen

- 1. Entfernen Sie alle Flussverbindungen zu anderen Systemmodulen:
 - Hinterspülpumpen-Einlassschlauch vom Peristaltikschlauch an der Hinterspülpumpe
 - Schläuche f
 ür das Waschen der Nadel (mit Beschriftung W und S) von den Degaser-Ports
 - Kapillare zum Autosampler vom Pumpenausgang
 - Lösungsmittelschläuche von den Pumpeneingängen

ACHTUNG

Flussverbindungen und Kapillaranschlüsse sind äußerst empfindlich gegen Verunreinigungen. Staub und Verschmutzungen können diese Anschlüsse verunreinigen.

Verschließen Sie Kapillaren stets mit Schutzkappen und offene Flussverbindungen mit Verschlussstopfen, um sie vor Verunreinigungen zu schützen.

- Verschließen Sie offene Flussverbindungen mit geeigneten Verschlussstopfen, zum Beispiel den Stopfen aus dem Zubehörkit.
- Kontrollieren Sie die Leckage-Wanne und den Drainage-Ablauf. Befindet sich Flüssigkeit darin, saugen Sie die Flüssigkeit mit einem Tuch auf.

7.20.2 Entfernen des Moduleinschubs der Pumpe



VORSICHT—Schweres und unhandliches Gerät

Die System-Module sind zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden an den System-Modulen zu vermeiden, beachten Sie folgende Hinweise:

- Um die Module zu greifen und zu transportieren, sind zwei Personen erforderlich, zum Beispiel, um sie anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn die Module in den Systemturm gehoben oder aus dem Systemturm entfernt werden.

Erforderliche Werkzeuge

Schraubendreher, Torx T20

Vorbereitungen

Siehe Vorbereiten für das Entfernen der Pumpe (> Seite 327).

Gehen Sie wie folgt vor

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



Abbildung 224: Unverlierbarkeitsschrauben am Moduleinschub der Pumpe

- 2. Drücken Sie alle Schläuche und Kapillaren, die sich im Schlauchkanal der Vanquish-Systemmodule befinden, in den Schlauchkanal hinein.
- 3. Greifen Sie den Moduleinschub an der Leckage-Wanne und ziehen Sie das Modul circa 10 cm aus dem Gehäuse heraus.

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.

7.20.3 Vorbereiten für das Entfernen des Autosamplers

Allgemeine Vorbereitungen

- 1. Nehmen Sie die Probenhalter und -gefäße aus dem Probenraum.
- 2. Führen Sie das Skript aus: **B05 Außerbetriebnahme (langfristig)** mit dem folgenden Parameter: **Zu entfernendes Modul = Autosampler**.
- 3. Soweit noch nicht geschehen, sichern Sie die Nadeleinheit (siehe Sichern der Nadeleinheit (▶ Seite 248)).

- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (> Seite 73).
- 5. Schieben Sie den Systemturm in eine geeignete Position.

Außerbetriebnahme

- Schalten Sie *alle* Module über ihren Hauptnetzschalter aus. Das Drücken des System-Ein- /Ausschalters reicht nicht aus, um die Geräte vollständig auszuschalten.
- 2. Ziehen Sie den Netzstecker.
- 3. Entfernen Sie alle Signalkabel zu anderen Geräten.

Flussverbindungen

- 1. Entfernen Sie alle Flussverbindungen zu anderen Systemmodulen und zu den Behältern:
 - Hinterspülschlauch und Einlassschlauch der Hinterspülpumpe. Verbinden Sie die Schläuche, die am Kopf der Dosiereinheit angeschlossen bleiben, mit einem Schlauchverbinder.
 - Nadel-Waschschläuche
 - Kapillare von der Pumpe
 - Kapillare zum Säulenthermostat, falls angeschlossen

ACHTUNG

Flussverbindungen und Kapillaranschlüsse sind äußerst empfindlich gegen Verunreinigungen. Staub und Verschmutzungen können diese Anschlüsse verunreinigen.

Verschließen Sie Kapillaren stets mit Schutzkappen und offene Flussverbindungen mit Verschlussstopfen, um sie vor Verunreinigungen zu schützen.

Reinigen

- 1. Wenn im Probenraum Kondensation oder verschüttete Proben vorhanden sind, reinigen und dekontaminieren Sie den Probenraum.
- 2. Kontrollieren Sie die Leckage-Wannen auf Flüssigkeiten hin. Tritt Flüssigkeit aus, saugen Sie die Flüssigkeit mit einem Tuch auf.

7.20.4 Entfernen des Moduleinschubs des Autosamplers



VORSICHT—Schweres und unhandliches Gerät

Die System-Module sind zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden an den System-Modulen zu vermeiden, beachten Sie folgende Hinweise:

- Um die Module zu greifen und zu transportieren, sind zwei Personen erforderlich, zum Beispiel, um sie anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn die Module in den Systemturm gehoben oder aus dem Systemturm entfernt werden.

Erforderliche Werkzeuge

Schraubendreher, Torx T20

Vorbereitungen

Siehe Vorbereiten für das Entfernen des Autosamplers (> Seite 329).

Entfernen des Moduleinschubs

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



Abbildung 225: Unverlierbarkeitsschrauben am Moduleinschub des Autosamplers

2. Drücken Sie alle Schläuche und Kapillaren, die sich im Schlauchkanal der Vanquish-Systemmodule befinden, in den Schlauchkanal hinein.

3. Greifen Sie den Moduleinschub am Mittelrahmen, wo sich das Türschloss befindet, und ziehen Sie das Modul circa 10 cm aus dem Gehäuse heraus.

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.

7.20.5 Vorbereiten für das Entfernen des Säulenthermostaten

Allgemeine Vorbereitungen

- Führen Sie das Skript aus: B05 Außerbetriebnahme (langfristig) mit dem folgenden Parameter: Zu entfernendes Modul = Säulenthermostat.
- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (> Seite 73).
- 3. Schieben Sie den Systemturm in eine geeignete Position.

Außerbetriebnahme

- 1. Schalten Sie *alle* Module über ihren Hauptnetzschalter aus. Das Drücken des System-Ein- /Ausschalters reicht nicht aus, um die Geräte vollständig auszuschalten.
- 2. Ziehen Sie den Netzstecker.
- 3. Entfernen Sie alle Signalkabel zu anderen Geräten.

Flussverbindungen

- 1. Entfernen Sie alle Flussverbindungen zu anderen Systemmodulen.
- 2. Versehen Sie offene Schaltventilöffnungen mit Verschlussfittingen.

ACHTUNG

Flussverbindungen und Kapillaranschlüsse sind äußerst empfindlich gegen Verunreinigungen. Staub und Verschmutzungen können diese Anschlüsse verunreinigen.

Verschließen Sie Kapillaren stets mit Schutzkappen und offene Flussverbindungen mit Verschlussstopfen, um sie vor Verunreinigungen zu schützen.

Optionales Zubehör

- Entfernen Sie die Säule gemäß der Anweisungen in Installieren von Säulen im Säulenthermostaten (▶ Seite 94) in umgekehrter Reihenfolge.
- 2. Entfernen Sie den Vorheizer, falls installiert (siehe Tauschen des passiven Vorheizers (▶ Seite 307)).
- Entfernen Sie die Schaltventile (siehe Tauschen des Vanquish-Schaltventils (▶ Seite 300) oder Tauschen des VICI-Schaltventils (▶ Seite 303)).

Reinigen

Kontrollieren Sie die Leckage-Wanne und den Drainage-Ablauf. Tritt Flüssigkeit aus, saugen Sie die Flüssigkeit mit einem Tuch auf.

7.20.6 Entfernen des Moduleinschubs des Säulenthermostaten

Erforderliche Werkzeuge

Schraubendreher, Torx T20

Vorbereitungen

Siehe Vorbereiten für das Entfernen des Säulenthermostaten (> Seite 332).

Entfernen des Moduleinschubs

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



Abbildung 226: Unverlierbarkeitsschrauben am Moduleinschub des Säulenthermostaten

- 2. Drücken Sie alle Schläuche und Kapillaren, die sich im Schlauchkanal der Vanquish-Systemmodule befinden, in den Schlauchkanal hinein.
- 3. Greifen Sie den Moduleinschub an der Leckage-Wanne und ziehen Sie das Modul circa 10 cm aus dem Gehäuse heraus.

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.

7.20.7 Zurückschicken des Moduleinschubs



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.

Vorbereitungen

Entfernen Sie den Moduleinschub, den Sie zurückschicken wollen, aus dem Gehäuse. Folgen Sie den Schritten in den jeweiligen Kapiteln:

Entfernen des Moduleinschubs der Pumpe (Seite 328)

Entfernen des Moduleinschubs des Autosamplers (Seite 331)

Entfernen des Moduleinschubs des Säulenthermostaten (Seite 333)

Gehen Sie wie folgt vor

- Bauen Sie den Moduleinschub in das Gehäuse des Ersatzteilgeräts ein. Folgen Sie den Schritten für den Einbau des Moduleinschubs in das Gehäuse und zum Anziehen der Unverlierbarkeitsschrauben in Installieren des Moduleinschubs (>>> Seite 336).
- 2. Wenden Sie sich an den Thermo Fisher Scientific-Kundendienst vor Ort, wenn Sie das Gerät für eine Reparatur an Thermo Fisher Scientific zurückschicken müssen.

7.20.8 Installieren des Moduleinschubs



VORSICHT—Schweres und unhandliches Gerät

Die System-Module sind zu schwer und zu unhandlich, um von nur einer Person sicher transportiert werden zu können. Um Verletzungen oder Schäden an den System-Modulen zu vermeiden, beachten Sie folgende Hinweise:

- Um die Module zu greifen und zu transportieren, sind zwei Personen erforderlich, zum Beispiel, um sie anzuheben oder zu verschieben.
- Mindestens zwei Personen sind insbesondere erforderlich, wenn die Module in den Systemturm gehoben oder aus dem Systemturm entfernt werden.

Erforderliche Werkzeuge

Schraubendreher, Torx T20

Vorbereitungen

1. 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 Neo-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 so weit ein, bis 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 227: Einsetzen des Moduleinschubs

- 4. Damit Sie den Moduleinschub leichter im Gehäuse befestigen können, ziehen Sie den Moduleinschub circa 5 mm Millimeter heraus.
- 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.
- Um das Modul anzuschließen, folgen Sie den Schritten in Installation (▶ Seite 75).
- 9. Lassen Sie die Seriennummer des Systems von einem Thermo Fisher Scientific Servicetechniker in das Systemmodul programmieren.

7.21 Vorbereiten des Systems für den Transport

ACHTUNG

Flussverbindungen und Kapillaranschlüsse sind äußerst empfindlich gegen Verunreinigungen. Staub und Verschmutzungen können diese Anschlüsse verunreinigen.

Verschließen Sie Kapillaren stets mit Schutzkappen und offene Flussverbindungen mit Verschlussstopfen, um sie vor Verunreinigungen zu schützen.

Allgemeine Vorbereitungen

- Führen Sie das Skript aus: B05 Außerbetriebnahme (langfristig) mit dem folgenden Parameter: Zu entfernendes Modul = Gesamtes System.
- 2. Soweit noch nicht geschehen, sichern Sie die Nadeleinheit (siehe Sichern der Nadeleinheit (▶ Seite 248)).
- Um besser an alle Systemkomponenten zu kommen, entsperren Sie den Systemsockel mit den Systemsockelschlüsseln. So können Sie den Systemsockel etwas auf dem Arbeitstisch verschieben. Siehe Entsperren des Systemsockels (>> Seite 73).
- 4. Schieben Sie den Systemturm in eine geeignete Position.
- 5. Nehmen Sie die Probenhalter und -gefäße aus dem Probenraum.

Lösungsmittel-System, Hinterspülung und Nadel-Waschsystem

TIPP Die Lösungsmittelschläuche müssen in der Regel nicht aus den Schlauchführungen entfernt werden (auch nicht, wenn der Moduleinschub getauscht wird). Wenn Sie jedoch die Lösungsmittelschläuche aus den Schlauchführungen in der Pumpe und allen Modulen oberhalb der Pumpe im Systemturm entfernen, achten Sie darauf, nicht an den anderen Schläuchen in den Führungen zu ziehen.

1. Leeren Sie alle Flüssigkeitsbehälter.

Außerbetriebnahme

 Schalten Sie *alle* Module über ihren Hauptnetzschalter aus. Das Drücken des System-Ein-/Ausschalters reicht nicht aus, um die Systemmodule vollständig auszuschalten.

	TIPP Wenn der Autosampler ausgeschaltet wird, öffnet sich die linke Tür des Autosamplers automatisch, um eine ordnungsgemäße Belüftung des Probenraums zu ermöglichen. Die Tür lässt sich nicht schließen, solange der Autosampler ausgeschaltet ist.		
	2.	Ziehen Sie alle Netzstecker.	
	3.	Entfernen Sie alle Signalkabel zu externen Geräten.	
Reinigen			
	1.	Wenn sich im Probenraum Kondenswasser gebildet hat oder dort Proben verschüttet wurden, reinigen und dekontaminieren Sie den Probenraum, bevor Sie fortfahren. Vergewissern Sie sich, dass der Autosampler ordnungsgemäß gereinigt und/oder dekontaminiert wurde.	
	2.	Kontrollieren Sie die Leckage-Wannen auf Flüssigkeiten hin. Tritt Flüssigkeit aus, saugen Sie die Flüssigkeit mit einem Tuch auf.	
Restliche Schritte			
	1.	Ziehen Sie den Ablaufschlauch aus dem Abfallbehälter.	
	2.	Bringen Sie die Tragegriffe an (siehe Anbringen der Tragegriffe (▶ Seite 340)), wenn nötig.	
	3.	Transportieren Sie das System.	
Inbetriebnehmen des Systems nach dem Transport			
	1.	Entsichern Sie die Nadeleinheit (siehe Entsichern der Nadeleinheit () Seite 249)).	
	2.	Schließen Sie die Kabel an (siehe Anschließen der Kabel (▶ Seite 107)).	
	3.	Schieben Sie den Systemturm zurück in Betriebsposition.	

- 4. Sperren Sie den Systemsockel mit den Systemsockelschlüsseln (siehe Sperren des Systemsockels (▶ Seite 118)).
- 5. Füllen Sie die Behälter (siehe entsprechende Abschnitte).
- 6. Stellen Sie die Flussverbindungen her (siehe Anschließen der Flussverbindungen (▶ Seite 119)).
- 7. Schalten Sie das System ein (siehe Einschalten des Systems (> Seite 149)).
- 8. Bereiten Sie das System für den Betrieb vor (siehe Vorbereiten des Systems für den ersten Betrieb (▶ Seite 138)).

7.22 Anbringen der Tragegriffe

Wann

Das System wird an einen anderen Ort transportiert.

Erforderliche Teile

4 Griffe

Verwenden Sie die Tragegriffe, die bei Auslieferung des Systems installiert waren und die 2 Bohrlöscher aufweisen.



Abbildung 228: Tragegriff für den Systemturm

• 8 Schrauben, die an den Griffen befestigt waren

Erforderliche Werkzeuge

Schraubendreher, Torx T30

Gehen Sie wie folgt vor



VORSICHT—Personenschäden und Schäden am System

Wenn Sie die Tragegriffe am Autosampler montieren und dann das System anheben, kann dies zu Personenschäden und zu Schäden am System führen.

Wenn Sie das gesamte System anheben und tragen müssen, befestigen Sie die Griffe nur an der Pumpe, um Personenschäden und Schäden am System zu vermeiden. 1. Schieben Sie die Tragegriffe von der Rückseite her in die rechte und linke Schiene.



Abbildung 229: Einschieben eines Tragegriffs in die rechte Schiene

2. Schieben Sie die Griffe so weit nach vorne, dass der Abstand zwischen der Vorderseite der Schiene und der Vorderseite des Griffes circa 6 cm beträgt.



Abbildung 230: Positionieren der vorderen Tragegriffe (obere Abbildung: Turm von links gesehen; untere Abbildung: Turm von rechts gesehen)

- 3. Ziehen Sie die Schrauben der Griffe an.
- 4. Schieben Sie die Tragegriffe von der Rückseite her in die rechte und linke Schiene.

- 5. Schieben Sie die Griffe so weit nach vorne, dass die Hinterseite der Griffe mit der Hinterseite der Schienen bündig abschließt.

Abbildung 231: Korrekte Position der hinteren Tragegriffe

6. Ziehen Sie die Schrauben der Griffe an.



VORSICHT—Verletzungsgefahr und Schäden am Modul

Die Tragegriffe können vom Modul herausrutschen, wenn die Befestigungsschrauben nicht korrekt angezogen sind. Außerdem kann ein Moduleinschub herausfallen, wenn er nicht korrekt befestigt ist. Dies kann zu Personenschäden und Schäden am Modul führen. Prüfen Sie Folgendes, bevor Sie das System anheben:

- Die Tragegriffe sind korrekt angeschraubt und können nicht aus dem Modul herausrutschen.
- Die vier Schrauben an allen Moduleinschüben sind festgezogen und kein Moduleinschub kann herausfallen (siehe Tauschen des Moduleinschubs (> Seite 327)).

8 Fehlersuche

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

8.1 Allgemeine Informationen zur Fehlersuche

An folgenden Dingen können Sie erkennen, dass während des Systembetriebs ein Problem aufgetreten ist.

Indikator	Beschreibung
Statusanzeigen	 Das System zeigt den Status folgender Komponenten an: System: Auf der Vanquish User Interface-Bildschirmseite oben. Siehe Systemstatus-Symbole (▶ Seite 155). Systemmodule Wenn die Firmware eines Systemmoduls ein Problem erkennt, erscheint eine Meldung in dem Benachrichtigungsfenster und die folgenden Statusanzeigen dieses Systemmoduls leuchten rot: LED-Statusleiste (siehe LED-Statusleisten (▶ Seite 153)) des
	 Moduls und die im Startfenster dargestellte LED-Statusleiste STATUS-LED (siehe Tastaturen () Seite 150))
Alarme bei Undichtigkeiten	Undichtigkeiten stellen ein potentielles Risiko für die Sicherheit und für das Chromatographie-Ergebnis dar. Wenn der Leaksensor auf eine Undichtigkeit anspricht, ertönt daher ein akustisches Signal, erscheint eine Meldung im Benachrichtigungsfenster und die Statusanzeige leuchtet rot. Folgen Sie den Anweisungen unter Testen auf Undichtigkeit () Seite 385).
Meldungen	Wenn die Firmware eines Systemmoduls ein Problem erkennt, erscheint eine Meldung in dem Benachrichtigungsfenster. Das Benachrichtigungsfenster speichert alle Warnungen und Fehler der Systemmodule und des System-Controllers seit dem letzten Booten des System-Controllers.
Analyseprobleme	Informationen zum Lösen von Analyseproblemen, wie zum Beispiel Peak- und Retentionszeitschwankungen, finden Sie in den entsprechenden Abschnitten unter System-Fehlersuche (> Seite 369).

In diesem Abschnitt finden Sie Informationen zu Meldungen und Störungen, die während des Systembetriebs auftreten können, und Sie finden Hilfe bei der Fehlersuche und Fehlerbehebung.

Zusätzliche Unterstützung

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. Halten Sie zur leichteren Identifizierung Ihres Gerätes bei der Kommunikation mit Thermo Fisher Scientific die Bestellnummer und die Seriennummer des Systems bereit. Diese Daten können Sie dem Typenschild entnehmen, das sich in der Schublade im Systemsockel befindet.

8.2 Meldungen

Wenn die Firmware eines Systemmoduls einen Fehlerzustand erkennt, erscheint eine Meldung im Vanquish User Interface und im Chromatographie-Datensystem. Die Tabellen in diesem Abschnitt führen die am häufigsten beobachteten Meldungen für jedes Systemmodul auf und leisten 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. Um die Ursache eines Problems weiter einzugrenzen, prüfen Sie, ob das Chromatographie-Datensystem zusätzliche Informationen zu einer bestimmten Meldung gibt.

Wenn Sie einen Code nicht finden können, prüfen Sie den Text der Meldung. Die beiden Meldungen "Unexpected module behavior" und "Module malfunction detected" können verschiedenen Codes zugewiesen sein.

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.

8.2.1 Übergreifende Meldungen

In diesem Abschnitt finden Sie Informationen zu den Meldungen, die für alle Systemmodule gelten.

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. 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.

Die Tabelle führt die am häufigsten beobachteten Meldungen für die Systemmodule auf.

Meldung und Code	Beschreibung und Abhilfe
Code 33 Leak detected – eliminate within approx. xx seconds.	Undichtigkeit erkannt - beheben Sie diese innerhalb von circa xx Sekunden.
	xx = Zeit in Sekunden, um die Undichtigkeit zu beheben, bevor die Sequenz abgebrochen wird
	Finden und beheben Sie die Ursache für die Undichtigkeit. Ist eine Undichtigkeit in der Pumpe nicht nach 3 Minuten behoben, stoppt der Pumpenfluss.
	 Unterziehen Sie alle Komponenten und Verbindungen im Flussweg einer optischen Pr
	 Ziehen Sie die undichte Verbindung oder Komponente nach oder tauschen Sie diese gegebenenfalls.
	3. Saugen Sie mit einem Tuch oder Papiertuch alle Flüssigkeit in der Leckage-Wanne und unter dem Leaksensor auf.
Code 34 Leak detected.	Undichtigkeit erkannt. Die Meldung erscheint, weil die Undichtigkeit, die zuvor mit Code 33 gemeldet wurde, nicht innerhalb von 3 Minuten behoben wurde. Aufgrund einer Undichtigkeit wurde die Pumpe gestoppt und/oder die laufende Sequenz abgebrochen. Finden und beheben Sie die Ursache für die Undichtigkeit (siehe Testen auf Undichtigkeit () Seite 385)).
Code 36 Download failed.	Download fehlgeschlagen.
	Der Firmware-Download wurde nicht erfolgreich abgeschlossen. 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 52 Module software incomplete. Download firmware (again).	Modul-Software unvollständig. Laden Sie die Firmware (nochmals) herunter.
	Die Firmware ist unvollständig, zum Beispiel, weil die Kommunikation zwischen dem Chromatographie-Datensystem und dem Modul während des Firmware-Downloads unterbrochen war. Wiederholen Sie den Download.
Code 79 [Device] was not found.	[Device] wurde nicht gefunden.
	Die Meldung kann für einen Drucksensor in einem Pumpenkopf oder für den Systemdrucksensor im Flowmeter erscheinen (siehe Informationen zur Identifikation der Ursache in Meldungen der Pumpe (▶ Seite 352)).
	 Pumpenkopf: Vergewissern Sie sich, dass das Drucksensorkabel des Pumpenkopfes, für den die Meldung erscheint, ordnungsgemäß angesteckt ist.
	• Flowmeter: Wenden Sie sich an den Kundendienst.

Meldung und Code	Beschreibung und Abhilfe
Code 84 Could not move X to its destination position. Code 85 Could not move X to its destination position.	X kann die Zielposition nicht anfahren. X = linkes Ventil oder rechtes Ventil Das entsprechende Schaltventil ist möglicherweise defekt. Informationen zum Austausch finden Sie unter Tauschen des Vanquish-Schaltventils (▶ Seite 300) oder Tauschen des VICI- Schaltventils (▶ Seite 303). Bleibt die Meldung bestehen, wenden Sie sich an den Kundendienst.
Code 89 Liquid leak sensor missing.	Leaksensor fehlt. Wenden Sie sich für Unterstützung an den Kundendienst von Thermo Fisher Scientific. Beachten Sie auch, dass das Modul dennoch betrieben werden kann, selbst wenn der Leaksensor defekt oder nicht installiert ist, indem der Leaksensor ausgeschaltet wird.
Code 90 Download firmware mismatch – invalid version.	Fehler bei Firmware-Download – ungültige Version. Sie haben versucht, eine inkompatible Firmware mit einer früheren Versionsnummer aufzuspielen als aktuell im Modul installiert ist. Das Aufspielen einer früheren Firmware-Version kann zu Funktionsverlusten oder Fehlfunktionen des Moduls führen. Falls erforderlich, wiederholen Sie die Installation mit einer Firmware-Version, die neuer ist als die aktuell im Modul installierte Version.
Code 91 Data changed during read.	 Datenänderung während des Lesevorgangs. Alle Module: Dies kann ein temporärer Fehler sein. Schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Modul wieder ein. Die Firmware ist möglicherweise defekt. Aktualisieren Sie die Firmware. Zusätzlich für die Pumpe: Wenn die Meldung für einen Drucksensor in einem Pumpenkopf oder für den Systemdrucksensor im Flowmeter erscheint (siehe Informationen zur Identifikation der Ursache in Meldungen der Pumpe (▶ Seite 352)): Pumpenkopf: Vergewissern Sie sich, dass das Drucksensorkabel des Teils, für das die Meldung erscheint, ordnungsgemäß angesteckt ist. Flowmeter: Wenden Sie sich an den Kundendienst.

Meldung und Code	Beschreibung und Abhilfe
Code 92 Data verification failed.	Datenverifikation fehlgeschlagen.
	Alle Module:
	 Dies kann ein temporärer Fehler sein. Schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Gerät wieder ein.
	 Die Firmware ist möglicherweise defekt. Aktualisieren Sie die Firmware.
	Zusätzlich für die Pumpe:
	Wenn die Meldung für einen Drucksensor in einem Pumpenkopf oder für den Systemdrucksensor im Flowmeter erscheint (siehe Informationen zur Identifikation der Ursache in Meldungen der Pumpe () Seite 352)):
	 Pumpenkopf: Vergewissern Sie sich, dass das Drucksensorkabel des Teils, für das die Meldung erscheint, ordnungsgemäß angesteckt ist.
	• Flowmeter: Wenden Sie sich an den Kundendienst.
Code 118 USB Buffer Overflow.	Überlauf des USB-Puffers.
	Es handelt sich um ein Software-Problem. Das Modul produziert Daten schneller als sie der Rechner, auf dem das Chromatographie-Datensystem läuft, verarbeiten kann.
	1. Trennen Sie das Modul im Chromatographie-Datensystem und verbinden Sie es erneut.
	 Wird das Problem damit nicht gelöst, aktualisieren Sie die Firmware oder die Version des Chromatographie-Datensystems.
	 Besteht das Problem weiterhin: Auch Software von Drittanbietern auf dem Rechner, zum Beispiel Virenscanner, oder eine schlechte Rechnerleistung, können das Problem verursachen. Wenden Sie sich an die IT-Abteilung vor Ort.
Code 120 System interlink request timed	Zeitüberschreitung der System-Interlink-Anforderung.
out.	Mit dem Modul kann keine Kommunikation aufgebaut werden. Das Modul hat nicht innerhalb der Zeit geantwortet.
	Für das Modul, für das die Meldung erscheint:
	1. Schalten Sie das Modul ein, wenn es noch nicht eingeschaltet ist.
	 Prüfen Sie die System-Interlink-Verbindungen zum Modul. Vergewissern Sie sich, dass alle System-Interlink-Kabel an beiden Enden angeschlossen sind.
	3. Bleibt die Meldung bestehen, tauschen Sie die System-Interlink- Kabel.

Meldung und Code	Beschreibung und Abhilfe
Code 126 Device lost system interlink connection.	 System-Interlink-Verbindung des Geräts wurde unterbrochen. Alle anderen Geräte am System-Interlink-Bus wurden ausgeschaltet oder ein Firmware-Download wurde auf allen Geräten durchgeführt. Prüfen Sie, ob alle anderen Module ausgeschaltet wurden oder ob auf allen anderen Modulen am Bus ein Firmware-Download durchgeführt wird. Warten Sie bis eines der Module am Bus wieder eingeschaltet wird oder bis der Firmware-Download beendet ist. Ein System-Interlink-Kabel ist möglicherweise vom Modul getrennt oder defekt. Prüfen Sie die System-Interlink-
	System-Interlink-Kabel an beiden Enden angeschlossen sind. 4. Bleibt die Meldung bestehen, tauschen Sie die System-Interlink-Kabel.
Code 128 System interlink data errors.	 System-Interlink-Datenfehler. Ein System-Interlink-Kabel ist möglicherweise vom Modul getrennt oder defekt. Prüfen Sie die System-Interlink-Verbindungen zum Modul. Vergewissern Sie sich, dass alle System-Interlink-Kabel an beiden Enden angeschlossen sind. Bleibt die Meldung bestehen, tauschen Sie die System-Interlink-
Code 129 Device X not found on system interlink.	 Gerät X am System-Interlink nicht gefunden. X = Seriennummer Das Modul mit der angegebenen Seriennummer ist möglicherweise nicht verbunden, ausgeschaltet oder die Firmware wird gerade aktualisiert. 1. Prüfen Sie die System-Interlink-Verbindungen zum Modul. 2. Bleibt die Meldung bestehen, prüfen Sie den Betriebszustand des Moduls.
Code 131 System interlink bus not operable.	 System-Interlink-Bus nicht betriebsbereit. Ein System-Interlink-Kabel ist möglicherweise vom Modul getrennt oder defekt. 1. Prüfen Sie die System-Interlink-Verbindungen zum Modul. Vergewissern Sie sich, dass alle System-Interlink-Kabel an beiden Enden angeschlossen sind. 2. Bleibt die Meldung bestehen, tauschen Sie die System-Interlink- Kabel.
Code 133 System Interlink access not supported.	Zugriff auf System-Interlink nicht unterstützt. Der Zugriff durch den System Controller ist nicht erlaubt. Sie können die Modul-Einstellungen nur über das Chromatographie- Datensystem ändern. Warten Sie zum Beispiel bis die Sequenz beendet ist und die LED- Leiste grün leuchtet. Dann wiederholen Sie den Vorgang über den System Controller.

Meldung und Code	Beschreibung und Abhilfe
Code 134 USB access by CDS not	USB-Zugriff durch CDS nicht unterstützt.
supported.	Der Zugriff durch das Chromatographie-Datensystem ist nicht erlaubt. Sie können die Modul-Einstellungen nur über den System Controller ändern.
	Warten Sie zum Beispiel bis das Skript beendet ist und die LED- Leiste grün leuchtet. Dann wiederholen Sie den Vorgang über das Chromatographie-Datensystem.
Code 136 Lock request rejected – already	Sperranforderung abgelehnt - bereits durch X gesperrt.
locked by X.	X = ID der sperrenden Komponente, mit Tasten-ID, USB-Adresse, die sich auf das Chromatographie-Datensystem bezieht oder System-Interlink-Adresse, die sich auf den System-Controller oder ein Modul bezieht
	Das Modul ist bereits durch eine andere Software (System- Controller oder Chromatographie-Datensystem) oder eine Taste gesperrt.
	Warten Sie, bis das Modul aus dem gesperrten Zustand freigegeben wird.
Code 137 Lock by X expired.	Sperre durch X ist abgelaufen.
	X = ID der sperrenden Komponente, mit Tasten-ID, USB-Adresse, die sich auf das Chromatographie-Datensystem bezieht oder System-Interlink-Adresse, die sich auf den System-Controller oder ein Modul bezieht
	Informieren Sie Thermo Fisher Scientific über das Auftreten. Es ist keine weitere Aktion erforderlich.
Code 145 Lock holder X lost.	Sperrende Komponente X verloren.
	X = ID der sperrenden Komponente, mit Tasten-ID, USB-Adresse, die sich auf das Chromatographie-Datensystem bezieht oder System-Interlink-Adresse, die sich auf den System-Controller oder ein Modul bezieht
	Die Sperre wurde automatisch freigegeben, da die sperrende Komponente vom System-Interlink verschwunden ist.
	1. Prüfen Sie, ob das sperrende Modul ausgeschaltet wurde.
	2. Prüfen Sie die System-Interlink-Verbindungen zum Modul.
Code 152 Assignment or command	Einstellung oder Befehl abgelehnt - Gerät gesperrt.
rejected – device locked.	Das Modul ist nicht bereit, Befehle oder Property-Einstellungen anzunehmen, da es noch von einer anderen Software gesperrt ist (System-Controller oder Chromatographie-Datensystem).
	Warten Sie, bis das Modul von der sperrenden Software freigegeben wird (System-Controller oder Chromatographie-Datensystem).

8.2.2 Meldungen der Pumpe

In diesem Abschnitt finden Sie Informationen zu den Meldungen mit den Code-Nummern 4000 bis 4999. Diese Meldungen sind spezifisch für die Pumpe.

Zusätzliche Identifikation der Ursache

Um die Ursache eines Problems weiter einzugrenzen, prüfen Sie die Meldung im Chromatographie-Datensystem auf zusätzliche Informationen:

Standardname im Datensystem	Beschreibung
PumpModule.Pump	Die Meldung bezieht sich auf das gesamte Pumpenmodul, einschließlich Flowmeter (mit Systemdrucksensor).
PumpModule.Pump Left Block (WP)	Die Meldung bezieht sich auf Folgendes:
	 • Linker Pumpenblock (A) • Komponenten im <i>linken</i> Pumpenblock, zum Beispiel im Pumpenkopf
	• Komponenten auf der <i>linken</i> Seite des <i>linken</i> Pumpenkopfs, zum Beispiel, auf den <i>linken</i> Drucksensor im Pumpenkopf oder den Arbeitskolben, der %A fördert.
PumpModule.Pump	Die Meldung bezieht sich auf Folgendes:
Left Block (EP)	• <i>Linker</i> Pumpenblock (A)
	 Komponenten im <i>linken</i> Pumpenblock, zum Beispiel im Pumpenkopf
	 Komponenten auf der rechten Seite des linken Pumpenkopfs, zum Beispiel, auf den rechten Drucksensor im Pumpenkopf oder den Ausgleichskolben, der %A fördert
PumpModule.Pump	Die Meldung bezieht sich auf Folgendes:
Right Block (WP)	• <i>Rechter</i> Pumpenblock (B)
	 Komponenten im <i>rechten</i> Pumpenblock, zum Beispiel im Pumpenkopf
	• Komponenten auf der <i>linken</i> Seite des <i>rechten</i> Pumpenkopfs, zum Beispiel, auf den <i>linken</i> Drucksensor im Pumpenkopf oder den Arbeitskolben, der %B fördert
PumpModule.Pump	Die Meldung bezieht sich auf Folgendes:
Right Block (EP)	• <i>Rechter</i> Pumpenblock (B)
	 Komponenten im <i>rechten</i> Pumpenblock, zum Beispiel im Pumpenkopf
	 Komponenten auf der rechten Seite des rechten Pumpenkopfs, zum Beispiel, auf den rechten Drucksensor im Pumpenkopf oder den Ausgleichskolben, der %B fördert

Meldungen der Pumpe

Die Tabelle führt die am häufigsten beobachteten Meldungen für die Pumpe auf.

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. 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 4098 Upper pressure limit	Obere Druckgrenze überschritten.
exceeded	Prüfen Sie den Flusspfad auf Durchlässigkeit. Führen Sie das Skript aus: D01 - Systemrückdruck testen .
Code 4099 Pressure fallen below lower	Der Druck liegt unter der unteren Druckgrenze.
limit.	Prüfen Sie die folgenden Ursachen:
	 Ein oder mehrere Lösungsmittelbehälter sind leer. Führen Sie das Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern (Nur aktualisieren = An).
	 Möglicherweise befinden sich Luftblasen in den Lösungsmittel- schläuchen. Kontrollieren Sie die Lösungsmittelfilter. Purgen Sie die Pumpe.
	 Bei dem eingestellten Fluss kann die Pumpe nicht genug Druck aufbauen, um die untere Druckgrenze zu erreichen. Passen Sie den unteren Grenzwert an oder erhöhen Sie den Fluss.
	 Möglicherweise gibt es eine Undichtigkeit im System. Führen Sie das Skript aus: E04 - Detaillierter Dichtigkeitstest (Pumpe).
	 Ein Kugelventil kann verschmutzt oder defekt sein. Reinigen oder tauschen Sie die Kugelventile. Siehe Warten der Pumpenkopf-Kugelventile (▶ Seite 224).
Code 4100 Purge pressure limit exceeded	Drucklimit für das Purgen überschritten
	Nach dem Start eines Purge-Zyklus hat sich ein Druck von mehr als 5 MPa (50 bar) aufgebaut.
	 Führen Sie das Skript aus: D01 - Systemrückdruck testen. Führen Sie die Abhilfen durch, die im Skript genannt werden.
	• Wenn die Meldung zusammen mit einer anderen Meldung zum Rückdruck des Flowmeter-Filters erscheint, (Code 4533), tauschen Sie den Filter aus (siehe Tauschen des Flowmeter- Filters () Seite 229)).
	 Das Purge-Ventil könnte defekt sein. Wenden Sie sich an den Kundendienst.

Meldung und Code	Beschreibung und Abhilfe
Code 4101 Not ready. Please run self- test.	Nicht bereit. Führen Sie einen Selbsttest durch. Die Pumpe ist nicht bereit, da der Selbsttest nicht erfolgreich war. Wiederholen Sie den Selbsttest. Führen Sie das Skript aus: B04 - Systemselbsttest .
Code 4112 Self-test failed because of fluctuating pressure.	Der Selbsttest ist aufgrund von Druckschwankungen fehl- geschlagen. Während des Selbsttests sind Druckschwankungen aufgetreten. Warten Sie, bis das System druckfrei ist oder machen Sie das System druckfrei und wiederholen Sie den Selbsttest. Führen Sie das Skript aus: B04 - Systemselbsttest .
Code 4113 Self-test failed because pressure is out of range.	Der Selbsttest ist fehlgeschlagen, da der Druck außerhalb des Bereichs liegt. Der Druck war während des Selbsttests nicht Null. Überprüfen Sie, dass die Kabel der Drucksensoren korrekt eingesteckt sind. Warten Sie, bis das System druckfrei ist oder machen Sie das System druckfrei und wiederholen Sie den Selbsttest. Führen Sie das Skript aus: B04 - Systemselbsttest . Wenn die Meldung erneut erscheint, ist gegebenenfalls der Systemdrucksensor defekt. Wenden Sie sich an den Kundendienst.
Code 4114 The device is busy. Please retry after 1 minute.	Das Gerät ist beschäftigt. Versuchen Sie es nach 1 Minute erneut. Die Pumpe ist nicht bereit, weil zum Beispiel ein Selbsttest oder eine Neukalibrierung läuft. Versuchen Sie es nach 1 Minute erneut. Wenn die Meldung erneut erscheint, warten Sie, bis das System druckfrei ist oder machen Sie das System druckfrei und schalten Sie die Pumpe aus und wieder ein.
Code 4119 Can't start pump while alarm is on.	Die Pumpe kann nicht gestartet werden, weil ein Alarm vorliegt. Es liegt ein Alarm vor, zum Beispiel, weil eine Undichtigkeit erkannt wurde. Sie können den Pumpenfluss erst dann erneut starten, wenn der Alarm nicht mehr vorliegt. Schalten Sie den Alarm stumm; drücken Sie dazu die Taste MUTE ALARM auf der Tastatur. Finden und beseitigen Sie die Ursache für den Alarm entsprechend den Angaben in der begleitenden Meldung.

Meldung und Code	Beschreibung und Abhilfe
Code 4120 The rear seal wash system has run out of wash solution.	Die Hinterspülung hat keine Waschlösung mehr.
	Diese Meldung warnt Sie nur, wenn die Tropfenzähler- funktionalität des Hinterspüldetektors eingeschaltet ist (Standardeinstellung).
	 Der Behälter für die Hinterspülflüssigkeit könnte leer sein. Stellen Sie neue Hinterspülflüssigkeit bereit und führen Sie das Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern (Hinterspülflüssigkeit hinten, Nur aktualisieren = An). Beachten Sie die Empfehlungen für die Hinterspülflüssigkeit in Füllen des Behälters für die Hinterspülflüssigkeit (> Seite 133).
	 Prüfen Sie die Hinterspülschläuche über den gesamten Flusspfad optisch auf Anzeichen von Blockade oder Undichtigkeit. Alle Schläuche müssen ordnungsgemäß angeschlossen und verlegt sein. Tauschen Sie gegebenenfalls die Hinterspülschläuche aus. Siehe Wartung der Hinterspülung (▶ Seite 279).
	 Prüfen Sie, ob Flüssigkeit an den Verbindungen der Hinterspülung oder unter den Pumpenköpfen austritt. Siehe Testen der Hinterspülung auf Undichtigkeit (> Seite 279).
	 Prüfen Sie den Peristaltikschlauch auf Anzeichen von Blockade oder Undichtigkeit. Tauschen Sie gegebenenfalls den Schlauch aus. Siehe Tauschen des Peristaltikschlauchs in der Hinterspülpumpe () Seite 289).
	 Überprüfen Sie, dass der Peristaltikschlauch korrekt in der Hinterspülpumpe eingelegt und der Hebel der Pumpe nicht blockiert ist.
	 Überprüfen Sie, dass der Hinterspülschlauch mit dem richtigen Port des Hinterspüldetektors verbunden ist (Port mit der Kennzeichnung in, siehe Bild in Testen der Hinterspülung auf Undichtigkeit (> Seite 279)).
	 Prüfen Sie die Elektroden des Hinterspüldetektors optisch auf Verschmutzung oder Beschädigung. Tauschen Sie gegebenenfalls den Hinterspüldetektor aus. Siehe Tauschen des Hinterspüldetektors () Seite 293).
Code 4121 Piston seal leakage has exceeded the recommended limit.	Die Undichtigkeit der Kolbendichtungen hat den empfohlenen Grenzwert überschritten.
	Diese Meldung warnt Sie nur, wenn die Tropfenzähler- funktionalität des Hinterspüldetektors eingeschaltet ist (Standardeinstellung).
	Testen Sie die Pumpe auf Dichtigkeit. Führen Sie das Skript aus: E04 - Detaillierter Dichtigkeitstest (Pumpe) . Wenn die Kolbendichtungen undicht sind, tauschen Sie die Dichtungen aus. Siehe Tauschen der Kolbendichtung (▶ Seite 208).

Meldung und Code	Beschreibung und Abhilfe
Code 4122 The rear seal leak sensor detects drops constantly.	Der Hinterspüldetektor detektiert permanent Tropfen.
	Diese Meldung warnt Sie nur, wenn die Tropfenzähler- funktionalität des Hinterspüldetektors eingeschaltet ist (Standardeinstellung).
	 Überprüfen Sie, dass die Flüssigkeit korrekt durch den Drainage- Ablauf rechts an der Leckage-Wanne in die Systemdrainage abläuft.
	 Prüfen Sie die Unterseite des Hinterspüldetektors: Entfernen Sie die Detektoreinheit der Hinterspülung aus der Pumpe und trennen Sie den Detektor vom Trichter. Trocknen Sie die Unterseite mit einem Tuch oder Papiertuch. Prüfen Sie die Detektorelektroden. Wenn diese sauber und unbeschädigt sind, setzen Sie Detektor und Trichter wieder zusammen und bauen Sie die Einheit wieder in die Pumpe ein. Siehe Tauschen des Hinterspüldetektors () Seite 293).
	 Wenn die Meldung erneut erscheint, könnte der Hinterspüldetektor defekt sein. Tauschen Sie den Hinterspüldetektor. Siehe Tauschen des Hinterspüldetektors (> Seite 293).
Code 4125 Degasser malfunction	Fehlfunktion des Degasers
	Die Überwachungsfunktion des Vakuumniveaus hat festgestellt, dass das Degaservakuum nicht ausreichend ist.
	Aktualisieren Sie die Firmware auf Version 2.41 oder höher.
Code 4127 The pump drive is still in undock position. Please execute a dock command.	Der Pumpenantrieb befindet sich noch in der Position Undock. Führen Sie einen Dock-Befehl aus.
	Diese Meldung warnt Sie während der Pumpenwartung (zum Beispiel, beim Tausch der Kolbendichtungen oder der Kolben), wenn Sie versuchen, die Pumpe zu starten, während sich der Pumpenkopf noch nicht in der korrekten Position für den Betrieb befindet.
	Bewegen Sie den Pumpenkopf zurück in die Position für den Betrieb: Die LEDs neben der Taste FLOW zeigen an, welcher Pumpenkopf sich in der Position für die Wartung befindet (LED blinkt grün). Vergewissern Sie sich, dass dieser Pumpenkopf gewählt ist (LED neben der Taste SELECT leuchtet grün) und drücken Sie die Taste DOCK . Wenn sich die Pumpe in der Position für den Betrieb befindet, blinkt die LED neben der Taste FLOW nicht mehr.
Code 4148 Can't perform this command while the flow is on.	Dieser Befehl kann nicht ausgeführt werden, während der Pumpenfluss angeschaltet ist.
	Sie haben versucht, einen Selbsttest durchzuführen, während der Fluss nicht null war. Schalten Sie den Fluss aus und wiederholen Sie den Befehl.

Meldung und Code	Beschreibung und Abhilfe
Code 4152 Pressure sensor malfunction. Check cable connection and retry.	Fehlfunktion des Drucksensors. Prüfen Sie die Kabelverbindung und versuchen Sie es erneut.
	Vergewissern Sie sich, dass das Kabel des Drucksensors, für den die Meldung erscheint, ordnungsgemäß angesteckt ist. Schalten Sie den Fluss wieder an.
	Wenn die Meldung erneut erscheint, ist gegebenenfalls der Drucksensor defekt:
	 Wenn die Meldung f ür den Drucksensor eines Pumpenkopfs erscheint, tauschen Sie den Pumpenkopf.
	 Wenn die Meldung f ür den Systemdrucksensor erscheint, wenden Sie sich an den Kundendienst.
Code 4158 Out of eluent X.	Kein Eluent X mehr vorhanden.
	Diese Meldung warnt Sie nur, wenn Sie den Lösungsmittel- verbrauch überwachen. In der Meldung ist das Lösungsmittel angegeben, auf das sich die Meldung bezieht. Stellen Sie neues Lösungsmittel zur 'Verfügung. Führen Sie das Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern .
Code 4161 Pressure recalibration	Abweichung des Drucks bei Neukalibrierung um xx bar.
deviates by xx bar.	Die Pumpe war während des Selbsttests nicht druckfrei. Warten Sie, bis die Pumpe druckfrei ist und führen Sie den Test erneut durch. Führen Sie das Skript aus: B04 - Systemselbsttest .
Code 4208 System pressure too high. Please relieve pressure and retry.	Zu hoher Systemdruck. Verringern Sie den Druck und versuchen Sie es erneut.
	Gehen Sie wie folgt vor:
	 Überprüfen Sie, dass die Kabel der Drucksensoren korrekt eingesteckt sind. Stecken Sie die Kabel gegebenenfalls neu an.
	 Prüfen Sie den Flusspfad auf Durchlässigkeit. Führen Sie das Skript aus: D01 - Systemrückdruck testen.
	3. Führen Sie einen Selbsttest durch. Führen Sie das Skript aus: B04 - Systemselbsttest .
	Wenn die Meldung erneut erscheint, ist gegebenenfalls der Systemdrucksensor defekt. Wenden Sie sich an den Kundendienst.
Code 4209 Could not build up enough pressure. Please check the fitting plug.	Es konnte nicht genug Druck aufgebaut werden. Prüfen Sie das Verschlussfitting.
	Die Pumpe konnte bei der Kalibrierung der Drucksensoren nicht genug Druck aufbauen. Dies kann folgende Ursachen haben:
	• Es könnte Luft im System sein. Purgen Sie die Pumpe. Führen Sie das Skript aus: C02 - Pumpe purgen .
	 Möglicherweise liegt eine Undichtigkeit vor. Führen Sie das Skript aus: E04 - Detaillierter Dichtigkeitstest (Pumpe).

Meldung und Code	Beschreibung und Abhilfe
Code 4213 Head pressure does not match system pressure. Execute the self- test or perform pressure recalibration.	 Der Druck im Kopf passt nicht zum Systemdruck. Führen Sie den Selbsttest durch oder starten Sie eine Neukalibrierung des Drucks. 1. Prüfen Sie den Flussweg vom Pumpenkopf zum Flowmeter- Einlass optisch auf Anzeichen von Blockierungen, insbesondere die Viper-Enden. Wenn Sie eine Blockierung vermuten, tauschen Sie die Kapillare. 2. Führen Sie einen Selbsttest durch. Führen Sie das Skript aus: B04 - Systemselbsttest. 3. Bleibt die Meldung bestehen, führen Sie eine Neukalibrierung des Drucks durch. Führen Sie das Skript aus: C05 - System- Drucksensoren abgleichen. 4. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Code 4220 Pump head not recognized. Please plug in the pressure sensor cable.	Pumpenkopf nicht erkannt. Stecken Sie das Kabel des Drucksensors an. 1. Stecken Sie das Drucksensorkabel vom Pumpenkopf am Port Pressure an. Wenn das Kabel angesteckt ist, überprüfen Sie es auf
	korrekten Sitz. 2. Wenn die Meldung weiterhin erscheint, könnte der Pumpenkopf defekt sein. Wenden Sie sich an den Kundendienst oder tauschen Sie den Pumpenkopf aus (siehe Tauschen des Pumpenkopfs (▶ Seite 195)).
Code 4231 Pressure sensor not recognized. Please plug in the sensor cable.	Drucksensor nicht erkannt. Stecken Sie das Sensorkabel an. Vergewissern Sie sich, dass das Kabel des Drucksensors, für den die Meldung erscheint, ordnungsgemäß angesteckt ist. Schalten Sie den Fluss wieder an.
	 Wenn die Meldung erneut erscheint, könnte der Drucksensor defekt sein: Wenn die Meldung für den Drucksensor eines Pumpenkopfs erscheint, tauschen Sie den Pumpenkopf (siehe Tauschen des Pumpenkopfs (> Seite 195)).
	 Wenn die Meldung f ür den Systemdrucksensor erscheint, k önnte der Systemdrucksensor defekt sein. Wenden Sie sich an den Kundendienst.
Code 4232 The identification number of the XX has changed.	Die Identifikationsnummer von XX hat sich geändert. XX ist der linke Pumpenkopf oder der rechte Pumpenkopf Die Meldung erscheint, wenn ein Pumpenkopf ausgetauscht wurde oder wenn der linke und der rechte Pumpenkopf vertauscht wurden, zum Beispiel versehentlich während Wartung/ Service. Die Meldung dient nur zur Information. Es sind keine Aktionen erforderlich.
Code 4234 The flow was stopped due to a request from another module.	Der Pumpenfluss wurde auf Anforderung eines anderen Moduls gestoppt. Der Pumpenfluss wurde aufgrund eines Problems bei einem anderen Modul gestoppt. Prüfen Sie die Meldungen von anderen Modulen.

Meldung und Code	Beschreibung und Abhilfe
Code 4238 Mainboard initialization failed.	 Fehler beim Initialisieren des Mainboards. Das Mainboard wurde nicht erfolgreich initialisiert. 1. Dies kann ein temporärer Fehler sein. Schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Gerät wieder ein. 2. Wenn die Meldung erneut erscheint, aktualisieren Sie die Firmware. 3. Wenn die Meldung weiterhin erscheint, könnte das Mainboard defekt sein. Wenden Sie sich an den Kundendienst.
Code 4503 The pressure exceeded the absolute limit.	Der Druck übersteigt den absoluten Grenzwert. Testen Sie den Flusspfad auf Durchlässigkeit. Führen Sie das Skript aus: D01 - Systemrückdruck testen .
Code 4504 NC drive initialization error.	Fehler bei der Initialisierung des NC-Antriebs. Wiederholen Sie den Selbsttest. Führen Sie das Skript aus: B04 - Systemselbsttest.
Code 4508 Piston drive home sensor remains dark.	Home-Sensor des Kolbenantriebs ist weiterhin dunkel. Wiederholen Sie den Selbsttest. Führen Sie das Skript aus: B04 - Systemselbsttest.
Code 4512 Cannot pressurize due to inlet valve failure or air in system.	 Druckbeaufschlagung nicht möglich, weil das Einlassventil defekt ist oder sich Luft im System befindet. Eine Kompressionsphase musste wiederholt werden, da der erwartete Druck nicht erreicht wurde. Es könnte sich Luft im System befinden oder das Einlasskugelventil muss gegebenenfalls gereinigt werden. Die Pumpe liefert jedoch den korrekten Fluss, da die folgende Kompressionsphase erfolgreich war. Wenn die Meldung selten ist und nicht zusammen mit anderen Meldungen erscheint, zum Beispiel Code 4532 oder Code 4538, können Sie sie tolerieren. Andernfalls: Überprüfen Sie den Füllstand im Lösungsmittelbehälter. Purgen Sie die Pumpe. Führen Sie das Skript aus: CO2 - Pumpe purgen. Wenn die Meldung nach dem Purgen erneut erscheint, reinigen Sie das Einlasskugelventil (siehe Reinigen der Kugelventile (Pumpenkopf) (▶ Seite 227)). Verwenden Sie nur Original-Lösungsmittelschläuche, die von Thermo Fisher Scientific für diese Pumpe autorisiert sind.
Code 4513 Home position deviation XX. Please execute self-test.	Abweichung der Home-Position XX. Führen Sie einen Selbsttest durch. XX gibt den Wert der Abweichung an. Vergewissern Sie sich, dass Sie die empfohlene Hinterspülflüssigkeit verwenden (siehe Füllen des Behälters für die Hinterspülflüssigkeit (> Seite 133)) und führen Sie einen Selbsttest durch. Führen Sie das Skript aus: B04 - Systemselbsttest .

Meldung und Code	Beschreibung und Abhilfe
Code 4514 The user defined calibration data of the flow sensor is invalid.	Die benutzerdefinierten Kalibrierdaten des Flusssenors sind ungültig.
	Die Kalibrierdaten für das ausgewählte Lösungsmittel sind unvollständig. Zum Beispiel, weil die Kalibrierprozedur unerwartet unterbrochen wurde.
	Wiederholen Sie die Kalibrierung für das ausgewählte Lösungs- mittel (siehe Hinzufügen von kundenspezifischen Lösungsmitteln zur Kalibriertabelle () Seite 168)).
Code 4517 Residual pressure too high for	Restdruck ist für das Fluss-Autozero zu hoch.
flow autozero.	Schalten Sie den Fluss und warten Sie, bis der Druck gefallen ist.
Code 4519 Calibration table full.	Kalibriertabelle voll.
	Löschen Sie ein Lösungsmittel aus der Kalibriertabelle der Lösungsmittel, damit Platz frei wird. Gehen Sie auf dem Skriptfenster zu Lösungsmittelkalibrierung. Führen Sie das Skript aus: M34 – Daten der Lösungsmittelkalibrierung löschen.
Code 4520 Flowmeter memory	Der Speicher des Flowmeters ist nicht verfügbar.
unavailable.	Überprüfen Sie, dass das Flowmeter korrekt eingebaut ist.
Code 4522 Unknown solvent.	Unbekanntes Lösungsmittel.
	Das von Ihnen gewählte Lösungsmittel ist nicht in der Kalibriertabelle der Lösungsmittel vorhanden. Mögliche Ursachen sind: Das Lösungsmittel wurde aus der Kalibriertabelle entfernt. Das Flowmeter wurde getauscht und ein Lösungsmittel (mit dem selben Namen) wurde nicht kalibriert und zu der Konfiguration des neuen Flowmeters hinzugefügt.
	Wählen Sie ein Lösungsmittel aus der Kalibriertabelle der Lösungsmittel oder fügen Sie das Lösungsmittel zur Kalibriertabelle hinzu (siehe Hinzufügen von kundenspezifischen Lösungsmitteln zur Kalibriertabelle (> Seite 168)).
Code 4523 Drift rate too high for flow	Drift-Rate ist für das Fluss-Autozero zu hoch.
autozero.	Der Fluss ist nicht stabil.
	1. Warten Sie, bis das System druckfrei ist.
	2. Purgen Sie die Pumpe, um Restluft zu entfernen.
	3. Wenn die Pumpe für mindestens 6 Stunden eingeschaltet war, wiederholen Sie den Autozero.
Code 4526 Command sequence error.	Fehler in der Befehlssequenz.
	Die Befehle wurden in einer falschen oder unerwarteten Reihenfolge ausgeführt. Prüfen Sie das Befehlsskript Ihrer Methode auf Fehler.
Code 4529 The pump capillary connections are setup for solvent calibration. Normal operation is not possible.	Die Kapillarverbindungen in der Pumpe sind für die Lösungsmittel- kalibrierung konfiguriert. Normalbetrieb ist nicht möglich.
	Ein Kalibrierversuch wurde unterbrochen. Verwenden Sie die Kalibrierskripte, um entweder die Kalibrierung abzuschließen oder um die ursprüngliche Kapillarkonfiguration der Pumpe entsprechend der Anwendung wiederherzustellen.
Meldung und Code	Beschreibung und Abhilfe
---	---
Code 4531 Cannot deliver flow at this pressure.	Der Fluss kann bei diesem Druck nicht geliefert werden. Die Pumpe kann die geforderte Flussrate bei dem sich daraus ergebenden Druck nicht liefern. Passen Sie die Anwendung an. Reduzieren Sie zum Beispiel die Flussrate. Siehe Abhängigkeit von verfügbarem Pumpendruck und Flussrate (▶ Seite 167).
Code 4532 Cannot regulate flow due to air in system or inlet valve failure.	 Fluss kann aufgrund von Luft im System oder defektem Einlassventil nicht geregelt werden. Eine oder mehrere Kompressionsphasen waren nicht erfolgreich. Die Pumpe konnte den korrekten Fluss nicht liefern. Es könnte sich Luft im System befinden oder das Einlasskugelventil muss gegebenenfalls gereinigt werden. Die Meldung erscheint zusammen mit Code 4512. Überprüfen Sie den Füllstand im Lösungsmittelbehälter. Purgen Sie die Pumpe. Führen Sie das Skript aus: C02 - Pumpe purgen. Wenn die Meldung nach dem Purgen erneut erscheint, reinigen Sie das Einlasskugelventil (siehe Reinigen der Kugelventile (Pumpenkopf) (▶ Seite 227)). Verwenden Sie nur Original-Lösungsmittelschläuche, die von
Code 4533 Elevated filter backpressure. Consider replacing the flowmeter filter soon.	Thermo Fisher Scientific für diese Pumpe autorisiert sind. Erhöhter Filterrückdruck. Tauschen Sie gegebenenfalls den Flowmeter-Filter bald. Die Meldung erscheint, bevor ein Problem auftritt. Die Meldung gibt an, dass es ein Problem mit den Lösungsmitteln geben könnte, zum Beispiel mit Algenwachstum. Tauschen Sie gegebenenfalls den Flowmeter-Filter (siehe Tauschen des Flowmeter-Filters (▶ Seite 229)). Beachten Sie die Empfehlungen unter Verwenden von Lösungsmitteln und Additiven (▶ Seite 145).
Code 4534 System pressure exceeded maximum pressure during purge.	 Systemdruck hat beim Purgen den maximalen Druck überschritten. 1. Entfernen Sie am Flowmeter-Ausgang die Kapillare zum Autosampler und versuchen Sie es erneut. 2. Ist wie folgt abhängig: Der Fehler tritt nicht mehr auf: Die Autosampler-Fluidik oder die Kapillare zum Autosampler könnten verstopft sein. Beheben Sie die Ursache für die Verstopfung und/oder tauschen Sie die Kapillare vom Flowmeter zum Autosampler aus. Der Fehler bleibt bestehen: Das Y-Stück oder die Widerstandskapillaren im Flowmeter könnten verstopft sein. Wenden Sie sich an den Kundendienst.
Code 4536 Solvent calibration failed.	 Lösungsmittelkalibrierung fehlgeschlagen. Möglicherweise sind Kapillarverbindungen oder Verschlussfittinge an den Pumpenköpfen oder am Flowmeter nicht korrekt befestigt oder undicht. 1. Ziehen Sie die Verbindungen, falls erforderlich, nach oder stellen Sie diese wieder her. 2. Wiederholen Sie die Kalibrierung.

Meldung und Code	Beschreibung und Abhilfe
Code 4538 Flow too low – check for	Fluss zu niedrig – prüfen Sie auf Undichtigkeiten.
leaks.	Die Pumpe kann die eingestellte Flussrate nicht liefern. Die Meldung kann alleine oder zusammen mit Code 4512 erscheinen.
	 Die Meldung erscheint zusammen mit Code 4512: Führen Sie die Abhilfemaßnahmen entsprechend der begleitenden Meldung aus. Wenn die Abhilfemaßnahmen nicht helfen, könnte eine Undichtigkeit im System vorliegen. Finden und beheben Sie die Ursache für die Undichtigkeit (siehe Testen auf Undichtigkeit (> Seite 385)).
	 Die Meldung erscheint alleine: Es könnte eine Undichtigkeit im System vorliegen. Finden und beheben Sie die Ursache für die Undichtigkeit (siehe Testen auf Undichtigkeit (▶ Seite 385)).
Code 4539 Flow too high.	Fluss zu hoch.
	Die Pumpe kann die Flussrate nicht auf den nominalen Fluss verringern.
	 Die Meldung erscheint, wenn die eingestellte Flussrate nahe Null ist (zum Beispiel, wenn der Pumpenfluss ausgeschaltet wird): Passen Sie die Druck- und Flusssensoren an. Führen Sie die Skripte aus: CO5 - System-Drucksensoren abgleichen und CO1 - Nullpunktabgleich des Pumpenflusssensors durchführen.
	 Die Meldung erscheint in einer anderen Situation: Wenden Sie sich an den Kundendienst.

8.2.3 Meldungen des Autosamplers

In diesem Abschnitt finden Sie Informationen zu den Meldungen mit den Code-Nummern 5000 bis 5999. Diese Meldungen sind spezifisch für den Autosampler.

Die Tabelle führt die am häufigsten beobachteten Meldungen für den Autosampler auf.

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. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Code 5002 Command rejected – module is in error state.	Befehl zurückgewiesen - Fehlerzustand im Modul aufgetreten. Ein interner Fehler ist aufgetreten. Führen Sie einen Selbsttest durch.
Code 5010 Horizontal needle drive blocked. Loosen transport lock completely.	Der horizontale Nadelantrieb ist blockiert. Lösen Sie die Transportsicherung vollständig. Lösen Sie die Transportsicherungsschraube oberhalb der Nadeleinheit. Siehe Entsichern der Nadeleinheit (* Seite 249). Prüfen Sie, ob der horizontale Nadelantrieb leicht verschoben werden kann. Wenn Hindernisse im Fahrweg des Nadelantriebs vorhanden sind, entfernen Sie diese.
Code 5011 Vertical needle drive blocked. Check for correct installation of sample loop.	Der vertikale Nadelantrieb ist blockiert. Prüfen Sie auf korrekten Einbau der Probenschleife. Prüfen Sie, ob die Probenschleife korrekt eingebaut ist. Folgen Sie den Schritten für den Einbau der Probenschleife in .
Code 5012 Selftest failed.	Selbsttest fehlgeschlagen. Der Autosampler-Selbsttest ist fehlgeschlagen. Prüfen Sie den Audit Trail auf zusätzliche Meldungen, die das Problem genauer spezifizieren. Wenden Sie sich für Unterstützung an den Kundendienst von Thermo Fisher Scientific.
Code 5013 Carousel blocked. Check for obstructions with racks or sample containers.	Der Drehteller ist blockiert. Prüfen Sie, ob Hindernisse durch Probenhalter oder Probengefäße bestehen. Drehen Sie den Drehteller und prüfen Sie, ob er sich leicht bewegen lässt. Wenn Hindernisse im Drehweg des Drehtellers vorhanden sind, entfernen Sie die Hindernisse.
Code 5014 Rack not found. Check for correct rack position in the carousel and sample sequence.	 Probenhalter nicht gefunden. Prüfen Sie, ob sich der Probenhalter in korrekter Position im Drehteller und in der Probensequenz befindet. Prüfen Sie, ob ein Probenhalter oder eine Mikrotiterplatte in diesem Segment im Drehteller eingesetzt sind. Prüfen Sie, dass Probenhalter bzw. Mikrotiterplatte korrekt eingesetzt sind (siehe) und korrigieren Sie die Positionierung, falls erforderlich.

Meldung und Code	Beschreibung und Abhilfe
Code 5015 Vial not found. Check for correct vial position in the rack and sample sequence.	Probenfläschchen nicht gefunden. Prüfen Sie, ob sich das Probenfläschchen in korrekter Position im Probenhalter und in der Probensequenz befindet.
	Prüfen Sie, ob sich ein Probenfläschchen oder eine Mikrotiterplatte an der angegebenen Position befindet. Stellen Sie ein Probengefäß an die angegebene Position (siehe). Prüfen Sie die festgelegten Probenpositionen in der Probensequenz. Legen Sie gegebenenfalls eine andere Position fest.
Code 5016 Needle seat not found. Install needle seat.	Nadelsitz nicht gefunden. Installieren Sie einen Nadelsitz. Vergewissern Sie sich, dass der Nadelsitz installiert ist. Installieren Sie den Nadelsitz und die Nadeleinheit oder tauschen Sie sie aus, falls erforderlich (siehe Tauschen des Nadelsitzes und der Nadeleinheit () Seite 251).
Code 5017 Needle hit an unexpected object due to incorrect sample rack or well plate configuration.	Die Nadel ist aufgrund einer inkorrekten Konfiguration von Probenhalter und Mikrotiterplatte unerwartet auf ein Objekt getroffen.
	Prüfen Sie, ob der Probenhalter bzw. die Mikrotiterplatte korrekt konfiguriert und die Probenposition korrekt in der Probensequenz definiert ist. Beachten Sie dabei die Anzahl der Positionen im konfigurierten Probenhalter bzw. in der konfigurierten Mikrotiterplatte sowie die jeweilige Plattenhöhe.
Code 5018 Vial pusher hit an object that is higher than expected. Check the rack and/or well plate configuration.	Der Niederhalter ist auf ein Objekt getroffen, das höher ist als erwartet. Prüfen Sie die Konfiguration von Probenhalter bzw. Mikrotiterplatte.
	Prüfen Sie, ob der richtige Probenhalter oder die richtige Mikrotiterplatte konfiguriert ist. Beachten Sie dabei die Höhe des Probenhalters bzw. der Mikrotiterplatte.
Code 5020 Septum too strong/tight for bottom detection. Check documentation	Septum zu stark/fest für Bodenerkennung. Der Dokumentation entnehmen Sie, welche Verbrauchsmaterialien geeignet sind.
for supported consumables.	Die Bodenerkennung funktioniert nicht mit dem aktuellen Septum. Verwenden Sie ein unterstütztes Septum oder schalten Sie die Bodenerkennung aus. Empfehlungen zu unterstützten Kappen und Verschlüssen von Probengefäßen finden Sie unter Empfehlungen zu Kappen und Verschlüssen von Probengefäßen (* Seite 166).
Code 5021 Insulation cover removed while temperature control is on. Replace cover or temperature control will shut	Die Isolierabdeckung wurde bei eingeschalteter Temperaturregelung entfernt. Tauschen Sie die Abdeckung aus oder die Temperatursteuerung schaltet in X.X Minuten ab.
down in X.X minutes.	Wobei X.X = die Dauer in Minuten
	Die (Probenschleifen-)Isolierabdeckung muss für Autosampler- Typen, die eine solche Abdeckung besitzen, montiert sein, damit die Probenraumthermostatisierung verwendet werden kann. Wenn die Abdeckung nicht ordnungsgemäß angebracht ist, kann die Thermostatisierung nicht eingeschaltet werden bzw. wird automatisch abgeschaltet, um den Innenraum vor hoher Feuchtigkeit und übermäßiger Kondensation zu schützen, wenn der Probenraum thermostatisiert wird.
	Wenn die Meldung erscheint, obwohl die Abdeckung montiert ist, prüfen Sie, dass der obere Teil der Abdeckung fest sitzt (im Bereich des Magneten). Sie können im Instrument Audit Trail nachsehen, ob die Abdeckung beim Abnehmen und wieder Aufsetzen korrekt erkannt wird.

Meldung und Code	Beschreibung und Abhilfe
Code 5022 Temperature control shut down due to missing loop cover.	 Abschaltung der Temperaturregelung wegen fehlender Probenschleifen-Isolierabdeckung. Die (Probenschleifen-)Isolierabdeckung muss für Autosampler- Typen, die eine solche Abdeckung besitzen, montiert sein, damit die Probenraumthermostatisierung verwendet werden kann. Wenn die Abdeckung nicht ordnungsgemäß angebracht ist, kann die Thermostatisierung nicht eingeschaltet werden bzw. wird automatisch abgeschaltet, um den Innenraum vor hoher Feuchtigkeit und übermäßiger Kondensation zu schützen, wenn der Probenraum thermostatisiert wird. Wenn die Meldung erscheint, obwohl die Abdeckung fest sitzt (im Bereich des Magneten). Sie können im Instrument Audit Trail nachsehen, ob die Abdeckung beim Abnehmen und wieder Aufsetzen korrekt erkannt wird.
Code 5023 Insulation Loop Cover Bracket removed while temperature control is on. Replace frame or temperature control will shut down in X.X minutes.	 Die Halterung der Probenschleifen-Isolierabdeckung wurde bei eingeschalteter Temperaturregelung entfernt. Setzen Sie die Halterung wieder ein oder die Temperaturregelung schaltet in X.X Minuten ab. Wobei X.X = die Dauer in Minuten Die Halterung für die Isolierabdeckung wurde entfernt während die Temperaturregelung eingeschaltet wurde. Die Isolierabdeckung muss montiert sein, um die Probenraumthermostatisierung zu verwenden. Bauen Sie die Halterung für die Isolierabdeckung wie in Tauschen des Nadelsitzes und der Nadeleinheit (▶ Seite 251) beschrieben wieder ein und installieren Sie die Isolierabdeckung wie in Einbauen der Isolierabdeckung (▶ Seite 99) beschrieben.
Code 5024 Insulation Loop Cover Bracket missing.	Halterung der Isolierabdeckung fehlt. Die Halterung für die Isolierabdeckung wurde entfernt während die Temperaturkontrolle eingeschaltet wurde. Bauen Sie die Halterung für die Isolierabdeckung wie in Tauschen des Nadelsitzes und der Nadeleinheit (> Seite 251) beschrieben wieder ein und installieren Sie die Isolierabdeckung wie in Einbauen der Isolierabdeckung (> Seite 99) beschrieben.
Code 5029 Carousel home sensor failure.	Home-Sensor des Drehtellers fehlgeschlagen. Wenn die Meldung erscheint, notieren Sie den Code der Meldung und schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Modul wieder ein. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Code 5030 Carousel motor failure or blockade.	Home-Sensor des Drehtellers fehlgeschlagen. Wenn die Meldung erscheint, notieren Sie den Code der Meldung und schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Modul wieder ein. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Code 5032 Both doors not closed while temperature control is on. Close both doors or temperature control will shut down in xx minutes.	Beide Türen bei eingeschalteter Temperaturregelung nicht geschlossen. Schließen Sie beide Türen oder die Temperaturregelung schaltet in xx Minuten ab. xx = Anzahl der Sekunden, bevor sich der Autosampler ausschaltet Halten Sie beide Autosamplertüren geschlossen, wenn die Temperaturregelung eingeschaltet ist.

Meldung und Code	Beschreibung und Abhilfe
Code 5033 Temperature control shut down due to open door.	Temperaturregelung wegen offener Tür abgeschaltet. Schließen Sie beide Autosamplertüren und schalten Sie die Temperaturregelung erneut ein.
Code 5035 Temperature control shut down due to xx fan failure.	Temperaturregelung wurde aufgrund eines xx Lüfterfehlers abgeschaltet. xx = Betroffener Lüfter Wenn die Meldung erscheint, notieren Sie den Code der Meldung und schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Modul wieder ein. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Code 5037 Backup memory error. Module not operational.	Fehler des Datensicherungsspeichers. Modul nicht betriebsbereit. Wenn die Meldung erscheint, notieren Sie den Code der Meldung und schalten Sie das Modul aus. Warten Sie 5 Sekunden und schalten Sie das Modul wieder ein. Wenden Sie sich an den Kundendienst, wenn die Meldung erneut erscheint.
Code 5039 Calibration process refused, pressure below 100 bar.	 Kalibriervorgang verweigert, Druck unter 100 bar. Drucksensorkalibrierung nicht möglich, da der Systemdruck zu gering ist. 1. Prüfen Sie das System auf Undichtigkeit, beschädigte Teile oder offene Verbindungen. 2. Purgen Sie den Autosampler: Führen Sie das Skript aus: C04 - Sampler purgen. 3. Prüfen Sie auf Undichtigkeiten oder Luft im System: Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler). 4. Wiederholen Sie die Drucksensorkalibrierung: Führen Sie das Skript aus: C05 - System-Drucksensoren abgleichen.
Code 5040 Model – Firmware Mismatch! This firmware only supports VN-A10 models! Download appropriate firmware!	Modell - Firmware stimmt nicht überein! Diese Firmware unterstützt nur VN-A10-Modelle! Laden Sie die richtige Firmware herunter! Sie haben versucht, die falsche Firmware herunterzuladen. Laden Sie eine Firmware-Datei für einen Vanquish Neo Autosampler herunter.
Code 5041 Abort error occurred. Command sequence aborted.	Abbruchfehler aufgetreten. Befehlssequenz abgebrochen. Prüfen Sie den Audit Trail auf zusätzliche Meldungen, die das Problem genauer spezifizieren.
Code 5042 Vanquish Neo sampler detected unknown inject valve bypass position. Set correct position and run self-test command.	Der Vanquish Neo-Sampler hat eine unbekannte Bypassposition des Injektionsventils erkannt. Stellen Sie die richtige Position ein und führen Sie den Selbsttestbefehl aus. Der Autosampler wurde während eines Skripts unterbrochen, wie z.B. dem Skript A06 - Fluidiksystem/Workflow ändern . Wiederholen Sie das Skript.
Code 5043 Vanquish Neo Pump not found on system interlink. Cannot receive pressure from pump SN xx.	Vanquish Neo Pumpe nicht im System-Interlink gefunden. Empfang des Drucks von Pumpe SN xx nicht möglich. Wobei xx = die Seriennummer der Pumpe Prüfen Sie, dass alle System-Interlink-Kabel am Vanquish-System korrekt angeschlossen sind.

Meldung und Code	Beschreibung und Abhilfe
Code 5044 Validation of vial bottom detection failed.	Die Validierung der Bodenerkennung ist fehlgeschlagen. Führen Sie den Befehl SelfTest aus oder schalten Sie den Autosampler aus und wieder ein, um die Validierung zu wiederholen.
Code 5045 Could not pressurize while delivering solvent. Check for air or leakages in system.	Druckbeaufschlagung während der Abgabe von Lösungsmittel nicht möglich. Prüfen Sie auf Luft oder Undichtigkeiten im System. Führen Sie folgendes Skript aus und folgen Sie den Anweisungen: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler).
Code 5099 The device is busy.	Das Gerät ist beschäftigt. Der Autosampler ist nicht bereit, weil zum Beispiel ein Selbsttest läuft oder der Autosampler noch einen vorherigen Befehl verarbeitet. Probieren Sie es erneut, wenn der Autosampler bereit ist.

8.2.4 Meldungen des Säulenthermostaten

In diesem Abschnitt finden Sie Informationen zu den Meldungen mit den Code-Nummern 2000 bis 2999. Diese Meldungen sind spezifisch für den Säulenthermostaten.

Die Tabelle führt die am häufigsten beobachteten Meldungen für den Säulenthermostaten auf.

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. 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 2095 Wrong right valve pod installed.	Falscher rechter Ventilkopf installiert. Prüfen Sie, dass der konfigurierte Ventilkopf und der installierte Ventilkopf übereinstimmen.
Code 2096 Wrong left valve pod installed.	Falscher linker Ventilkopf installiert. Prüfen Sie, dass der konfigurierte Ventilkopf und der installierte Ventilkopf übereinstimmen.
Code 2097 Right valve position error.	 Positionsfehler des rechten Ventils. 1. Prüfen Sie, dass der konfigurierte Ventilkopf und der installierte Ventilkopf übereinstimmen. 2. Prüfen Sie, dass der Ventilkopf korrekt installiert ist. 3. Starten Sie die Funktion "Catch Valve", indem Sie die Taste VALVE RIGHT auf der Tastatur gedrückt halten.
Code 2098 Left valve position error.	 Positionsfehler des linken Ventils. 1. Prüfen Sie, dass der konfigurierte Ventilkopf und der installierte Ventilkopf übereinstimmen. 2. Prüfen Sie, dass der Ventilkopf korrekt installiert ist. 3. Starten Sie die Funktion "Catch Valve", indem Sie die Taste VALVE LEFT auf der Tastatur gedrückt halten.

8.3 System-Fehlersuche

In diesem Abschnitt finden Sie einen Leitfaden zur Fehlersuche für den Betrieb des Vanquish Neo-Systems.

Er gibt Informationen zu Fehlerbildern und Ursachen, die direkt mit den LC-Systemmodulen zusammenhängen. Entnehmen Sie Informationen zur Fehlersuche in der Chromatographie im Allgemeinen und bei speziellen Applikationen der Fachliteratur.

Das Vanquish User Interface bietet einige Skripte zur Suche von Fehlern im System.

Probleme bei der Detektion

Problem	Skript
 Hoher Säulen-Rückdruck Überdruckabschaltung Schlechte Präzision der Peak- flächen Kontinuierlich steigender Druck Kein Fluss oder die Pumpe gibt eine Meldung aus, nach der sie den Fluss nicht regeln kann 	D01 - Systemrückdruck testen
 Niedriger Säulendruck Retentionszeitschwankungen Schlechte Präzision der Peak- flächen Kein Fluss oder die Pumpe gibt eine Meldung aus, nach der sie den Fluss nicht regeln kann 	D02 - Systemdichtigkeit testen (was getestet werden soll = System)

Probleme während der Analyse

Problem	Skript
Die Injektion oder Sequenz wurde abgebrochen und das automatisch gestartete B03 - System reinigen-Skript schlägt fehl	Führen Sie das Skript aus: B03 - System reinigen .

Sonstige Probleme

Problem	Skript
Fehlerzustand: Eine oder mehrere LED-Leisten leuchten rot und das Vanquish User Interface zeigt an, dass sich das System in einem Fehlerzustand befindet.	 Führen Sie das Skript aus: B04 - Systemselbsttest. Prüfen Sie, ob der Systemstatus ok ist (Symbol ist grün). Falls nicht, prüfen Sie das Benachrichtigungsfenster auf Meldungen und beheben Sie das Problem (siehe Meldungen () Seite 346)).
Systemeinstellungen passen nicht zueinander	Um beim System oder bei einzelnen Systemmodulen die werksseitigen Standardwerte wiederherzustellen, führen Sie das Skript aus: E02 – Auf System- Werkseinstellungen zurücksetzen.
Die Systemeinstellungen konnten nicht abgerufen und das System konnte nicht korrekt initialisiert werden	Führen Sie das Skript mit den für Ihre Konfiguration und ihren Workflow geeigneten Parameter durch: E01 - Systemeinrichtung starten.
Das Ändern des Fluidiksystems oder des Workflows (A06 - Fluidiksystem/Workflow ändern) wurde nicht erfolgreich abgeschlossen und nur wenige Skripte stehen zur Verfügung	

Suchen Sie in den folgenden Abschnitten 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.

8.3.1 Peaks

Keine Peaks

Mögliche Ursache	Abhilfe
Vorheizer wird in Nano- oder Kapillar-LC- Anwendungen ver- wendet	Entfernen Sie den Vorheizer (siehe Tauschen des passiven Vorheizers () Seite 307)).
Zwischen Säule und Detektor oder zwischen Autosampler und Säule wurde die falsche Kapillare ver- wendet	Verwenden Sie die für Ihre jeweilige Anwendung korrekten Kapillare (um den Workflow und die Flussverbindungen zu wechseln, führen Sie das Skript aus: A06 - Fluidiksystem/Workflow ändern oder entnehmen Sie diese Informationen aus dem <i>Flow Scheme</i>).
Luftblasen in der Probenschleife, Nadel oder Dosiereinheit	Purgen Sie den Autosampler.
Probenfläschchen leer	Überprüfen Sie den Füllstand im Probenfläschchen. Wenn der Füllstand zu niedrig ist, ersetzen Sie das Probenfläschchen durch ein aufgefülltes Probenfläschchen.
Lösungsmittel im Lösungsmittelbehälter entspricht nicht den Lösungsmitteleinstel- lungen	Verwenden Sie das Lösungsmittel, dass Sie eingestellt haben oder passen Sie die Einstellung an das Lösungsmittel im Lösungsmittelbehälter an.

Peakabweichung

Mögliche Ursache	Abhilfe
Temperatur des Säulen- thermostaten nicht stabil	 Prüfen Sie die Temperatur des Säulenthermostaten. Warten Sie, bis sich die Temperatur stabilisiert hat.
Nachdem der Lösungs- mitteltyp geändert wurde, wurde das vor- herige Lösungsmittel nicht vollständig aus dem Autosampler oder System herausgespült	 Wenn das Skript (B01 - Flüssigkeiten/Lösungsmittel ändern) das vorherige Lösungsmittel nicht vollständig herausgespült hat, prüfen Sie, ob der korrekte Parameter gesetzt wurde: Nur aktualisieren = Aus. Führen Sie das Skript erneut aus: B01 - Flüssigkeiten/Lösungsmittel ändern mit dem Parameter: Nur aktualisieren = Aus.

8.3.1.1 Peakform

Peak-Tailing

Mögliche Ursache	Abhilfe
Extrasäulenvolumen zu groß	 Verwenden Sie kurze Kapillaren mit geeignetem Innendurchmesser. Verwenden Sie die für Ihre Anwendung geeigneten Viper-Kapillare (Informationen hierzu finden Sie im Skript: A06 - Fluidiksystem/Workflow ändern).
Ungeeignete Kapillarver- bindungen	 Prüfen Sie, dass die Fittinge korrekt sitzen. Lösen und ziehen Sie die Viper-Fittinge an (siehe Anschließen von NanoViper- Kapillaren (▶ Seite 121)). Verwenden Sie die für Ihre Anwendung geeigneten Viper-Kapillare (Informationen hierzu finden Sie im Skript: A06 - Fluidiksystem/Workflow ändern). Tauschen Sie die Kapillare.
Totvolumina in den Ka- pillarverbindungen	 Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler). Überprüfen Sie die Kapillare und tauschen Sie sie aus, falls erforderlich.

Peakverbreiterung

Mögliche Ursache	Abhilfe
Extrasäulenvolumen zu groß	 Wenn frühere eluierende Peaks breiter sind als spätere, prüfen Sie zum Beispiel den Kapillar-Innendurchmesser und die Kapillarlänge.
	 Verwenden Sie eine Kapillare mit kleinerem Volumen f ür den Anschluss zwischen dem S
Temperaturgradient in der Säule	• Verwenden Sie bei Flussraten von > 50 μ L/min, insbesondere bei Säulen mit einem Innendurchmesser von 1 mm, in einem Flussbereich von 50-100 μ L/min einen Vorheizer.
	 Wenn bereits ein Vorheizer installiert ist, pr üfen Sie, ob dieser korrekt installiert ist. Falls erforderlich, installieren Sie den Vorheizer erneut (siehe Tauschen des passiven Vorheizers (Seite 307)).
Kapillare teilweise ver-	1. Führen Sie das Skript aus: D01 - Systemrückdruck testen.
stopft	2. Tauschen Sie die Kapillaren aus.
Probenschleife teil- weise verstopft	Führen Sie das Skript aus: D01 - Systemrückdruck testen .
Totvolumina in den Ka- pillarverbindungen	1. Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler).
	2. Befestigen Sie die Kapillarverbindungen korrekt.
	3. Ziehen Sie die Kapillarverbindungen nach.
	4. Tauschen Sie die Kapillare.

Peakverbreiterung, hohe Totzeit

Mögliche Ursache	Abhilfe
Nachsäulenvolumen hoch	 Verwenden Sie eine Kapillare mit den folgenden geeigneten Parametern: Innendurchmesser Länge Prüfen Sie die Kapillare. Wenn Sie defekt ist, tauschen Sie die Kapillare aus.
Nadelsitzfilter verstopft	 Führen Sie das Skript aus: D01 - Systemrückdruck testen. Tauschen Sie gegebenenfalls den Nadelsitz aus.

Dreieckpeaks

Mögliche Ursache	Abhilfe
Säule ist überladen (Konzentration der Pro- be zu hoch)	Verdünnen Sie die Probe.

8.3.1.2 Störpeaks, Negative Peaks und Spikes

Störpeaks

Mögliche Ursache	Abhilfe
Kontamination (typischerweise Injektor oder Säule)	Spülen Sie das System inklusive der Säule mit einem geeigneten Lösungsmittel (siehe Spülen bei hartnäckiger Verschmutzung (▶ Seite 189)).
Die verwendeten Eluenten sind alt oder verunreinigt oder die Qualität ist nicht aus- reichend	 Verwenden Sie frische und geeignete Lösungsmittel, zum Beispiel in MS-Qualität. Ersetzen Sie die Lösungsmittelbehälter durch Behälter mit frischen und geeigneten Lösungsmitteln (in MS-Qualität)(führen Sie das Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern).

Negative Peaks

Mögliche Ursache	Abhilfe
Lösungsmittel der Probe und der mobilen Phase unterscheiden sich in der Zusammensetzung	Lösen Sie die Probe in der mobilen Phase.
Luft injiziert/Proben- fläschchen leer	Prüfen Sie, ob das Probenfläschchen leer ist. Ersetzen Sie das Probenfläschchen durch ein aufgefülltes Probenfläschchen.

Spikes

Mögliche Ursache	Abhilfe
Es treten elektrische Störungen von anderen Systemmodulen auf	 Betreiben Sie keine Geräte in der Nähe, die starke elektrische Felder oder Magnetfelder erzeugen. Schließen Sie keine Geräte, die Stromschwankungen verursachen könnten, an dasselbe Stromnetz an, das auch das System mit Strom versorgt.
Säulentemperatur deutlich über dem Siedepunkt der mobilen Phase	Verhindern Sie bei der Methodenentwicklung, dass die Säulentemperatur deutlich über dem Siedepunkt der mobilen Phase liegt.
Lösungsmittel nicht ge- nügend entgast	Entgasen Sie die Lösungsmittel in einem Ultraschallbad.

8.3.1.3 Präzision der Peakflächen

Schlechte Präzision der Peakflächen/Peakflächen zu klein

Mögliche Ursache	Abhilfe
Luftblasen in der Probenschleife, Nadel oder Dosiereinheit	Purgen Sie den Autosampler.
Probenfläschchen leer	Überprüfen Sie den Füllstand im Probenfläschchen. Wenn der Füllstand zu niedrig ist, ersetzen Sie das Probenfläschchen durch ein aufgefülltes Probenfläschchen.

Kontamination oder Verschleppung

Mögliche Ursache	Abhilfe
Kontamination im Sys- tem	Spülen Sie das System mit einem geeigneten Lösungsmittel. Wenn Sie im Pumpenkanal A Wasser als Lösungsmittel verwenden, spülen Sie das System mit 80% Acetonitril.
Verschleppung des Autosamplers durch	1. Prüfen Sie, ob einer der Nadel-Waschschläuche (mit der Beschriftung SWP und WWP) gequetscht ist.
verunreinigte Nadel	 Prüfen Sie, ob einer der Nadel-Waschbehälter leer ist. Ist einer der Behälter leer, tauschen Sie ihn durch Ausführen des folgenden Skripts aus: B01 - Flüssigkeiten/ Lösungsmittel ändern. Verwenden Sie eine geeignete Nadel-Waschflüssigkeit (siehe Füllen der Behälter für die Nadel-Waschflüssigkeit (> Seite 130)).
	 Pr üfen Sie, ob das Waschen der Nadel im Chromatographie-Datensystem deaktiviert ist. Wenn ja, aktivieren Sie das Waschen der Nadel.
	 Um die Nadel und den Nadelsitz zu waschen, f ühren Sie das Skript aus: B03 - System reinigen.
	 Tritt die Verschleppung weiterhin auf, tauschen Sie die Nadeleinheit und den Nadelsitz aus (siehe Tauschen des Nadelsitzes und der Nadeleinheit (▶ Seite 251)).
Verschleppung des Autosamplers durch verschlissenen oder be- schädigten Nadelsitz	Tauschen Sie den Nadelsitz aus (siehe Tauschen des Nadelsitzes und der Nadeleinheit (▶ Seite 251)).

Analyse ist instabil oder nicht reproduzierbar

Mögliche Ursache	Abhilfe
Instabile Umgebungs- bedingungen	 Sorgen Sie f ür gleichm
	Thermostatisieren Sie die Säule.
	 Wenn die Probe gekühlt werden kann, kühlen Sie sie im Autosampler.
	• Vermeiden Sie Zugluft.
Gradient nicht reproduzierbar	 Prüfen Sie den %B-Wert. Wenn er nicht zwischen 1 und 99 liegt, korrigieren Sie ihn.
	 Das System ist nicht genügend äquilibriert. Lassen Sie es 5 Stunden lang äquilibrieren.
	• Purgen Sie die Pumpe.

Undichtigkeit

Mögliche Ursache	Abhilfe
Kapillarverbindungen nicht korrekt befestigt oder undicht	 Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = System). Überprüfen Sie die Kapillare und tauschen Sie sie aus, falls erforderlich.

Probenprobleme

Mögliche Ursache	Abhilfe
Probe ist instabil und zerfällt	 Kühlen Sie die Probe im Autosampler. Verwenden Sie eine neue Probe. Ändern Sie die Umgebungsbedingungen.

Injektionsvolumen variiert

Mögliche Ursache	Abhilfe
Autosampler saugt Luft an	Es ist zu wenig Probe im Probenfläschchen vorhanden oder die Nadelhöhe ist nicht korrekt eingestellt.
	 Überprüfen Sie den Füllstand im Probenfläschchen. Wenn der Füllstand zu niedrig ist, ersetzen Sie das Probenfläschchen durch ein aufgefülltes Probenfläschchen.
	 Um zu pr üfen, ob die Nadel verstopft ist, f ühren Sie das Skript aus: D01 - Systemr ückdruck testen. Folgen Sie den Anweisungen.
Luft im Flussweg des Autosamplers	Purgen Sie den Autosampler.
Gasgehalt der Probe zu hoch oder gesättigt	Entgasen Sie die Probe, wenn möglich.
Injektionsnadel ver- stopft oder Nadelspitze verformt	Führen Sie das Skript aus: D01 - Systemrückdruck testen . Folgen Sie den Anweisungen.
Kolbendichtungen in Autosampler-Dosierein- heit undicht	Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler) . Folgen Sie den Anweisungen.

Mögliche Ursache	Abhilfe
Injektionsventil oder andere Autosampler- Komponenten undicht	Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler) . Folgen Sie den Anweisungen.
Nadelsitzverschleiß	Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = Sampler) . Folgen Sie den Anweisungen.

8.3.2 Druck

Mögliche Ursache	Abhilfe
Das System ist mit Partikeln verstopft	Es stellt sich heraus, dass häufig Proben und ungenügende Lösungsmittelreinheit für Systemverstopfungen verantwortlich sind. Für Nano-LC- und Kapillar-LC- Anwendungen mit ihren noch kleineren Innendurchmessern werden eine noch höhere Lösungsmittel- und Probenreinheit benötigt.
	• Nano-LC- und Kapillar-LC-Anwendungen: Um Probenverunreinigungen zu vermeiden, beachten Sie Folgendes:
	 ♦ Verwenden Sie die Probenfläschchen mit partikelfreien Septen aus dem Zubehörkit oder die in Empfehlungen zu Kappen und Verschlüssen von Probengefäßen (▶ Seite 166) aufgeführten Probenfläschchen.
	 Reduzieren Sie die Anzahl von Replikaten aus einem Probenfläschchen mit Kappe auf <10.
	 Waschen Sie die Nadel und/oder intensivieren Sie das Nadelwaschen im Chromatographie-Datensystem.
	 Führen Sie das Skript aus: D01 - Systemrückdruck testen. Folgen Sie den Anweisungen.
Lösungsmittel ist ver- unreinigt	Verwenden Sie ein frisches Lösungsmittel.

Hoher Rückdruck

Mögliche Ursache	Abhilfe
Kapillaren im System verstopft oder geknickt	Führen Sie das Skript aus: D01 - Systemrückdruck testen . Folgen Sie den Anweisungen.
Säule verunreinigt oder verstopft	 Führen Sie das Skript aus: D01 - Systemrückdruck testen. Folgen Sie den Anweisungen. Wenn der Test keine Verstopfung feststellen kann, ist die Säule oder die Kapillare zum nachgeschalteten Detektionsgerät verstopft. Tauschen Sie die Säule oder die Kapillare zum nachgeschalteten Gerät.
Säulenschaltventil ver- stopft	 Prüfen Sie das Ventil auf Durchlässigkeit. Tauschen Sie das Ventil, falls erforderlich (siehe Tauschen des Vanquish- Schaltventils (> Seite 300) oder Tauschen des VICI-Schaltventils (> Seite 303)).
Vorheizer verstopft	Tauschen Sie den Vorheizer (siehe Tauschen des passiven Vorheizers (> Seite 307)).

Mögliche Ursache	Abhilfe
Kontamination im Sys- tem	Spülen Sie das System durch Purgen der Pumpe und des Autosamplers mit einem geeigneten Lösungsmittel.
Nadelsitz verstopft	Wenn das Rückdrucktest-Skript die Meldung ausgibt, dass im Nadelsitz ein hoher Rückdruck besteht, führen Sie das Skript aus: C21 - Nadeleinheit und Nadelsitz reinigen oder tauschen mit dem Parameter: Nadelsitz reinigen.
Lösungsmittel im Lösungsmittelbehälter entspricht nicht den Lösungsmitteleinstel- lungen	Verwenden Sie das Lösungsmittel, dass Sie eingestellt haben oder passen Sie die Einstellung an das Lösungsmittel im Lösungsmittelbehälter an.

Niedriger Rückdruck

Mögliche Ursache	Abhilfe
Undichtigkeit im System	Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = System) . Folgen Sie den Anweisungen.
Luft im Autosampler	Führen Sie das Skript aus: CO4 - Sampler purgen.
Lösungsmittel im Lösungsmittelbehälter entspricht nicht den Lösungsmitteleinstel- lungen	Verwenden Sie das Lösungsmittel, dass Sie eingestellt haben oder passen Sie die Einstellung an das Lösungsmittel im Lösungsmittelbehälter an.

8.3.3 Retentionszeitschwankungen

Abnehmende Retentionszeiten

Mögliche Ursache	Abhilfe
Ungeeignete Zusam-	 Prüfen Sie das Alter des vorgemischten Lösungsmittels. Purgen Sie die Pumpe. Wenn Sie das System für den Betrieb vorbereiten, lassen Sie es für 5-24 Stunden
mensetzung des	äquilibrieren. Führen Sie das Skript aus: C01 - Nullpunktabgleich des Pumpenflusssensors
Lösungsmittels	durchführen.

Ansteigende Retentionszeiten

Mögliche Ursache	Abhilfe
Ungeeignete Zusam- mensetzung des Lösungsmittels	 Prüfen Sie das Alter des vorgemischten Lösungsmittels. Purgen Sie die Pumpe. Wenn Sie das System für den Betrieb vorbereiten, lassen Sie es für 5-24 Stunden äquilibrieren. Führen Sie das Skript aus: C01 - Nullpunktabgleich des Pumpenflusssensors
	durchführen.

Mögliche Ursache	Abhilfe
Abnehmende Flussrate	1. Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = System). Folgen Sie den Anweisungen.
	2. Überprüfen Sie die eingestellte Flussrate.

Streuende Retentionszeiten

Mögliche Ursache	Abhilfe
Druckschwankungen aus Pumpe	Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = System).
Luft im Autosampler	Führen Sie das Skript aus: CO4 - Sampler purgen mit dem Parameter: Waschen des Nadelinneren, schwache Flüssigkeit.

Retentionszeitschwankungen über einen längeren Zeitraum hinweg

Mögliche Ursache	Abhilfe
Nach der Umstellung auf andere Lösungs- mittel wurde das vor- herige Lösungsmittel nicht vollständig aus der Pumpe heraus- gespült.	Purgen Sie die Pumpe.
Säulen-, Proben- oder Lösungsmittel- temperatur instabil	Warten Sie, bis die Säulen-, Proben- oder Lösungsmitteltemperatur vollständig äquilibriert ist.

Retentionszeit außerhalb des erwarteten Bereichs

Mögliche Ursache	Abhilfe
Die thermischen Flusssensoren haben eventuell eine kleine Drift oder eine Ver- schiebung.	Führen Sie das Skript aus: C01 - Nullpunktabgleich des Pumpenflusssensors durchführen.
Die vorgemischten Lösungsmittel sind zu alt.	 Verwenden Sie einen frischen Eluenten. Verwenden Sie frische und geeignete Lösungsmittel (zum Beispiel in MS-Qualität). Purgen Sie die Pumpe.
Vorheizer wird in Nano- LC-Anwendungen ver- wendet	Entfernen Sie den Vorheizer (siehe Tauschen des passiven Vorheizers () Seite 307)).
Nachsäulenvolumen falsch	Prüfen Sie das Volumen der Kapillare, die an den OUT-Port der Säule angeschlossen ist.
Falsche Probenschleife eingestellt	Prüfen Sie, ob die Einstellungen für die Probenschleife zu der installierten Probenschleife passen und korrigieren Sie die Einstellungen, falls erforderlich.

Mögliche Ursache	Abhilfe
Lösungsmittel im Lösungsmittelbehälter entspricht nicht den Lösungsmitteleinstel- lungen	Verwenden Sie das Lösungsmittel, dass Sie eingestellt haben oder passen Sie die Einstellung an das Lösungsmittel im Lösungsmittelbehälter an.

8.3.4 Basislinie

Starke Basisliniendrift

Mögliche Ursache	Abhilfe
System nicht ausrei- chend äquilibriert	 Spülen Sie das System, bis ein stabiles Gleichgewicht erreicht ist. In der Regel ist dies nach 5 - 10 Säulenvolumina der Fall.
	 Wird das Problem nicht durch das Sp ülen des Systems gelöst, purgen Sie die Pumpe und lassen Sie das System mindestens f ür 5 Stunden äquilibrieren.
Instabile Umgebungs- bedingungen	Sorgen Sie für gleichmäßige Temperatur und Luftfeuchtigkeit.
Die Absorption des Eluenten verändert sich, wenn der Gradient gefahren wird	Absorbierende Additive können das Absorptionsspektrum abhängig vom Lösungsmittel verändern. Erwägen Sie variierende Additivkonzentrationen, um die Drift auszugleichen.
Autosampler verunrei- nigt	Führen Sie das Skript aus: B03 - System reinigen .
Säule verunreinigt	Spülen oder ersetzen Sie die Säule.
Rückstände des alten Lösungsmittels im Kopf der Dosiereinheit oder in der Probenschleife	Um die Dosiereinheit oder die Probenschleife zu spülen, führen Sie das Skript aus: C04 - Sampler purgen mit dem Parameter: Waschen des Nadelinneren, starke Flüssigkeit . Um Lösungsmittelbehälter auszutauschen, führen Sie immer dieses Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern .
Die verwendeten Lö- sungsmittel sind verun- reinigt oder nicht ho- mogen	 Stellen Sie sicher, dass die Lösungsmittel in den Behältern vor der Analyse vollständig durchmischt sind. In wässrigen Lösungsmitteln können sich Mikroorganismen vermehren, welche das Lösungsmittel verunreinigen. Um den Lösungsmittelbehälter gegen einen Lösungsmittelbehälter mit frischem Lösungsmittel auszutauschen, führen Sie das Skript aus: B01 - Flüssigkeiten/Lösungsmittel ändern. Folgen Sie den Anweisungen.

Unregelmäßige Schwankungen der Basislinie, starkes Rauschen

Mögliche Ursache	Abhilfe
Luft im System	Purgen Sie die Pumpe und den Autosampler.
Das Lösungsmittel ist verunreinigt oder die Qualität ist nicht ausreichend	 Verwenden Sie einen frischen Eluenten. Verwenden Sie geeignete Lösungsmittel (MS-Qualität). Um auf ein anderes Lösungsmittel umzustellen, führen Sie die Skripte aus: A01 - Typen der Pumpenlösungsmittel einstellen und dann B01 - Flüssigkeiten/ Lösungsmittel ändern.

Periodische Schwankungen der Basislinie, Pulsation

Mögliche Ursache	Abhilfe
Druckschwankungen aus Pumpe	 Purgen Sie die Pumpe. Führen Sie das Skript aus: D01 - Systemrückdruck testen.
Luft im System	Purgen Sie die Pumpe.

8.3.5 Längere Ladezeit der Trap-Säule

Mögliche Ursache	Abhilfe
Die Trap-Säule ist veraltet oder verstopft	1. Führen Sie das Skript aus: D01 - Systemrückdruck testen mit dem Parameter: Nur Trap-Säulen-Widerstand testen.
	 Notieren Sie den Widerstandswert und vergleichen Sie ihn mit einem früheren Ergebnis.
	 Wenn der Widerstand zu hoch f ür Ihre Anwendung ist, f ühren Sie den gesamten R ückdrucktest aus. F ühren Sie das Skript aus: D01 - Systemr ückdruck testen.
	4. Überprüfen Sie das Testergebnis und tauschen Sie gegebenenfalls die verstopfte Komponente.
	 Wenn Sie eine Komponente getauscht haben, führen Sie den Rückdrucktest erneut aus: D01 - Systemrückdruck testen.
Nadelsitz verstopft	Siehe Druck (> Seite 376), Unterabschnitt Hoher Rückdruck.

8.3.6 Temperaturregelung

8.3.6.1 Temperaturregelungsprobleme beim Säulenthermostaten

Säulenthermostat-Temperatur ändert sich über längere Zeit nicht, obwohl der Sollwert noch nicht erreicht ist

Abhilfe
 Sorgen Sie für eine niedrigere Umgebungstemperatur (z.B. durch Klimatisieren des Raums). Wählen Sie einen höheren Temperatur-Sollwert.
Schalten Sie die Temperaturregelung ein.
 Öffnen Sie die Tür des Säulenthermostaten und warten Sie, bis sich die Lüfter nicht mehr drehen. Prüfen Sie, ob die Lüftungsöffnungen des Säulenthermostaten verdeckt sind. Entfernen Sie sämtliche Gegenstände vor den Lüftungsöffnungen

Mögliche Ursache	Abhilfe
Umgebungsluft dringt in Säulenraum ein	Eine unzureichende Abdichtung des Säulenraums kann die Heizleistung des Säulenthermostaten beeinträchtigen. 1. Prüfen Sie, ob die Gerätetüren korrekt schließen.
	 Prüfen Sie, ob die Gerätetüren richtig abdichten. Prüfen Sie, ob die Türdichtung beschädigt ist. Wenn die Türdichtung beschädigt ist, tauschen Sie die Gerätetür (siehe Tauschen der Türen (* Seite 325)).
Säulenthermostat defekt	Wenden Sie sich an den Kundendienst.

8.3.6.2 Temperaturregelungsprobleme beim Autosampler

Autosampleı	[.] erreicht	die Sollt	temperatur	nicht
-------------	-----------------------	-----------	------------	-------

Mögliche Ursache	Abhilfe
Umgebungstemperatur oder -luftfeuchtigkeit zu hoch	 Stellen Sie sicher, dass die Isolierabdeckung eingebaut ist. Sorgen Sie für eine niedrigere Umgebungstemperatur (z.B. durch Lüften) bzw. Luftfeuchtigkeit.

8.3.7 Stromversorgung und Kommunikation (Systemmodule)

Die in diesem Abschnitt beschriebenen Probleme und Abhilfen beziehen sich auf die Systemmodule (zum Beispiel Systemsockel, Pumpe, Autosampler und Säulenthermostat (wenn vorhanden)). Informationen zu Stromversorgungs- und Kommunikationsproblemen des System Controllers, Displays und der Benutzeroberfläche finden Sie unter Probleme bei System Controller, Display und der Benutzeroberfläche (\triangleright Seite 383).

Das assamts System hat keinen Strom (alle LED Leisten sind aus (dunk	
1 α β α β α β	- 11
	211
D a b	- 1 / /

Mögliche Ursache	Abhilfe
System-Ein-/Aus- schalter ausgeschaltet	Schalten Sie das System über den unteren-Ein-/Ausschalter vorne links am Systemsockel ein.

Status-LED am System-Controller ist aus (dunkel)

Mögliche Ursache	Abhilfe
Ein-/Ausschalter des System Controllers ist ausgeschaltet	Schalten Sie den System Controller über den oberen Ein-/Ausschalter vorne links am Systemsockel ein.
Hauptnetzschalter des System Controllers ist ausgeschaltet	Schalten Sie den System-Controller über den Hauptnetzschalter ein.
Gerät nicht am Stromnetz angeschlossen	Schließen Sie das Netzkabel an.
Fehler in der Elektronik	Wenden Sie sich an den Kundendienst.

Systemmodul hat keinen Strom (LED-Leiste aus (dunkel))

Mögliche Ursache	Abhilfe
Gerät ausgeschaltet	Schalten Sie das betreffende Gerät über den Hauptnetzschalter ein.
Gerät nicht am Stromnetz angeschlossen	Schließen Sie das Netzkabel an.
Sicherung durchgebrannt	Tauschen Sie die Sicherung aus (siehe Tauschen der Sicherungen () Seite 324)).
Fehler in der Elektronik	Wenden Sie sich an den Kundendienst.

Ein oder mehrere Geräte lassen sich nicht über den System-Ein-/Ausschalter am Systemsockel ausschalten

Mögliche Ursache	Abhilfe
System-Interlink-Kabel fehlen oder sind defekt	Prüfen Sie die System-Interlink-Verbindungen zwischen dem Systemsockel und den Systemgeräten, siehe Anschließen der System-Interlink-Kabel (Seite 113).

Modul lässt sich nicht vom Chromatographie-Datensystem steuern

Mögliche Ursache	Abhilfe
Keine Verbindung zwischen Gerät und Rechner	 Prüfen Sie die USB-Verbindungen vom System zum Rechner, siehe Anschließen der USB-Kabel (> Seite 109). Wenn das Gerät über den USB-Hub eines anderen Geräts angeschlossen ist, stellen Sie sicher, dass das Gerät mit dem Hub eingeschaltet ist. Überprüfen Sie den USB-Port am Rechner. Es wird USB-Standard 2.0 benötigt.

8.3.8 Probleme bei System Controller, Display und der Benutzeroberfläche

Fehlerbild	Mögliche Ursache	Abhilfen
Bild auf dem Display flackert	Das Display ist nicht korrekt verbunden	Prüfen Sie, ob das Display-Kabel fingerfest am Display Data -Port des System Controllers angezogen ist.
Display funktioniert nicht	Display nicht an den System Controller angeschlossen	 Prüfen Sie, ob das Display am System Controller angeschlossen ist. Wenn das nicht der Fall ist, gehen Sie zum nächsten Schritt über. Schalten Sie den System Controller aus. Schließen Sie das Display an den System Controller an (siehe Anschließen der Display-Kabel (> Seite 111)). Schalten Sie den System Controller ein.
System Controller kann nicht über seinen Ein-/Aus- schalter ausge- schaltet werden	Das kurze VSC-Kabel ist nicht verbunden	Stellen Sie sicher, dass das kurze VSC-Kabel den VSC- Port des Systemsockels mit dem VSC-Port des System Controllers verbindet.
System Controller ist eingeschaltet, aber die Status-LED am System- Controller ist aus (dunkel)		
System Controller kann nicht eingeschaltet werden, obwohl der Ein-/Aus- schalter gedrückt wurde		
Keine Skripte verfügbar	-	Schalten Sie den System Controller aus, warten Sie 5 Sekunden und schalten Sie ihn wieder ein.
Kein Systemmodul im Vanquish User Interface angezeigt - oder -	Systemturm nicht an den System-Interlink-Port des System Controllers angeschlossen	Schließen Sie den Systemturm an den System-Interlink- Port auf der kurzen Seite des System Controllers an (siehe Anschließen der System-Interlink-Kabel () Seite 113)).
Skripte fehlen im Vanquish User Interface	Systemturm nicht an den korrekten System-Interlink- Port des System Controllers angeschlossen	Schließen Sie den Systemturm an den System-Interlink- Port auf der kurzen Seite des System Controllers an (siehe Anschließen der System-Interlink-Kabel () Seite 113)).
Einzelnes Systemmodul nicht im Vanquish User Interface angezeigt - oder - Skripte fehlen im Vanquish User Interface	Einzelnes Systemmodul ist nicht an die System- Interlink-Schnittstelle ange- schlossen	 Prüfen Sie, ob das fehlende Systemmodul an die System-Interlink-Schnittstelle des Systems angeschlossen ist. Ist dies nicht der Fall, integrieren Sie es in den System-Interlink-Bus des Systemturms (siehe Anschließen der System-Interlink-Kabel (> Seite 113)). Prüfen Sie, ob das Systemmodul über den Hauptnetzschalter eingeschaltet ist. Tauschen Sie das System-Interlink-Kabel.

Fehlerbild	Mögliche Ursache	Abhilfen
	Firmware-Version des Moduls und/oder des System Controllers nicht korrekt oder Firmware- Versionen passen nicht zueinander	Installieren Sie die korrekten und passenden Firmwareversionen auf Ihrem Modul und/oder System Controller. Wenden Sie sich für Unterstützung an den Kundendienst von Thermo Fisher Scientific.
Seriennummer des Systems im Vanquish User Interface ist Nicht definiert	Nach dem Ersetzen eines Moduleinschubs wurde die Seriennummer nicht gesetzt.	Lassen Sie die Seriennummer des Systems von einem Thermo Fisher Scientific Servicetechniker in das Systemmodul programmieren.

8.4 Testen auf Undichtigkeit

Testen Sie regelmäßig, ob bei folgenden Flüssigkeitssystemen Undichtigkeiten im System vorhanden sind:

Flüssigkeit	Folgen Sie den Anweisungen in
Lösungsmittel	Diesem Abschnitt
Hinterspülflüssigkeit	Testen der Hinterspülung auf Undichtigkeit (▶ Seite 279)
Lösungsmittel und Nadel- Waschflüssigkeit von den Autosampler- Ventilen	Prüfen der Autosampler-Ventile auf Dichtigkeit (▶ Seite 296)
Kondenswasser aus dem Autosampler	Testen des Drainagepumpen-Schlauchs (▶ Seite 265)

Wann

- Regelmäßig (als Bestandteil des Wartungszeitplans):
 - Nach dem Tauschen der Lösungsmittelbehälter
 - Nach dem Wechseln des Workflows
- Ein Leaksensor hat eine Undichtigkeit erkannt.
- Bei chromatographischen Problemen:
 - Peaks sind zu klein
 - Retentionszeitschwankungen

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 177).

Gehen Sie wie folgt vor

1. Ist wie folgt abhängig:

Wann	Vorgehen
Test wird regelmäßig durchgeführt (als Bestandteil des Wartungszeitplans)	Führen Sie das Skript aus: D02 - Systemdichtigkeit testen (was getestet werden soll = System).
Test wird durchgeführt, weil	1. Führen Sie das Skript aus: D02 -
• ein Leaksensor eine Undichtigkeit	Systemdichtigkeit testen (was getestet
erkannt hat.	werden soll = System).
 chromatographische Probleme	 Führen Sie das Skript aus: E04 -
beobachtet werden.	Detaillierter Dichtigkeitstest (Pumpe).

- 2. Prüfen Sie das Testergebnis. Wenn das Skript eine Undichtigkeit erkannt hat, beheben Sie sie, indem Sie den Anweisungen auf dem Bildschirm folgen.
- 3. Um zu prüfen, ob die Undichtigkeit komplett behoben wurde, führen Sie das Skript erneut aus und prüfen Sie anschließend das Ergebnis.

9 Spezifikationen

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

9.1 Technische Spezifikationen

9.1.1 Umgebungsbedingungen

Die Umgebungsbedingungen 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
Überspannungskategorie	Ш

9.1.2 Leistungsaufnahme

Die Leistungsaufnahme ist wie folgt spezifiziert:

Produkt	Leistungsaufnahme
Pumpe	100 – 240 V AC, ± 10 %; 50/60 Hz, max. 525 W / 550 VA
Autosampler	100 – 240 V AC, ± 10 %; 50/60 Hz; max. 525 W / 550 VA
Säulenthermostat	100 – 240 V AC, ± 10 %; 50/60 Hz; max. 525 W / 550 VA
System Controller	100 – 240 VAC, ± 10 %; 50/60 Hz, max. 50 W / 150 VA

9.1.3 Abmessungen

Die Abmessungen des Systems und der Systemmodule sind wie folgt spezifiziert:

Produkt	Abmessungen (Höhe x Breite x Tiefe)	
System	Basiskonfiguration ¹ : 74 x 45 x 65 cm (Display am Autosampler befestigt; Position vor dem Solvent Rack)	Basiskonfiguration und Säulenthermostat: 82 x 45 x 65 cm (Display am Autosampler befestigt; Position vor dem Säulenthermostat)
Pumpe	19,2 x 42 x 62 cm	
Autosampler	29 x 42 x 62 cm	
Säulenthermostat	Außenmaße	15,9 x 42 x 62 cm
	Säulenraum	7,8 x 34,3 x 3,7 cm
¹ : Zur Basiskonfiguration gehört der Systemsockel mit System-Controller, Pumpe, Autosampler und das Display		

9.1.4 Emissionsschalldruckpegel

Die Emissionsschalldruckpegel sind wie folgt spezifiziert:

Produkt	Emissionsschalldruckpegel
System	< 75 dB(A), typisch < 55 dB(A)
Pumpe	< 70 dB(A), typisch < 50 dB(A)
Autosampler	< 70 dB(A), typisch < 50 dB(A)
Säulenthermostat	< 42 dB(A), typisch < 40 dB(A)

9.1.5 Gewicht

Das Gewicht des Systems und der Systemmodule ist wie folgt spezifiziert:

Produkt	Gewicht	
System	Basiskonfiguration ¹ : Circa 66 kg	Basiskonfiguration und Säulenthermostat: Circa 79 kg
Pumpe	24 kg	
Autosampler	26 kg	
Säulenthermostat	Ohne Ventile: 10,7 kg; mit 2 Vanquish-Ventilen: 11,6 kg; mit 2 VICI-Ventilen: 13,1 kg	
¹ : Zur Basiskonfiguration gehört der Systemsockel mit System Controller, Pumpe, Autosampler und das Display		

9.2 Leistungsspezifikationen

9.2.1 Leistungsspezifikationen des Systems

Art	Spezifikation
Kommunikation	Am Systemsockel befinden sich folgende Ports:
	1 System Interlink-Port (RJ45-8 Stecker)
	1 VSC-Port (RJ-12 Stecker)
	Informationen zu den Ports der Systemmodule finden Sie in den Abschnitten zur Leistungsspezifikation der einzelnen Systemmodule.
Steuerung	Vanquish User Interface
	Tastatur mit Funktionstasten zur Ausführung bestimmter Funktionen direkt am Systemmodul.
	Das System kann mit einem Chromatographie-Datensystem gesteuert werden. Weitere Fragen beantwortet Ihnen gern die Thermo Fisher Scientific-Vertriebsorganisation.
Materialien im Flussweg	Informationen zu den Materialien im Flussweg finden Sie in den Abschnitten zur Leistungsspezifikation der Pumpe, des Autosamplers und des Säulenthermostaten.
Informationen zu Lösungsmitteln und Additiven	Siehe Verwenden von Lösungsmitteln und Additiven () Seite 145).
Biokompatibilität	Ja
Sicherheitsmerkmale	Leakerkennung und sichere Leakbehandlung, Überdruckabschaltung
Good Laboratory Practice (GLP)	Predictive Performance-Funktionen zur Planung von Wartungsarbeiten basierend auf den tatsächlichen Betriebs- und Nutzungsbedingungen der Systemmodule.
	Alle Systemparameter werden im Benachrichtigungsfenster protokolliert.

Die Leistungsdaten des Systems sind wie folgt spezifiziert:

9.2.2 Leistungsspezifikationen der Pumpe

Art	Spezifikation
Funktionsprinzip	Serielle Zweikolbenpumpe mit unabhängigen Kolbenantrieben
Kompensation der Kompression	Vollautomatisch und unabhängig von der Zusammensetzung der mobilen Phase
Flussbereich (einstellbar)	1 nL/min – 100 μL/min in Schritten von 1 nL/min
Flussgenauigkeit	Typisch: ± 3% Bedingungen: 0,1 - 100 μL/min jedes der 4 vorkalibrierten Lösungsmittel, gemessen bei 20 MPa Systemdruck
Abweichung zwischen den Modulen	< 5% bei Flussraten > 300 nL/min – 10 μL/min Bedingungen: Gemessen durch Retentionszeitdifferenzen des Proteinverdau-Peptids TGQAPGFSYTDANK Cytochrom C und 30 min- Gradient.
Retentionszeitpräzision	≤ 0,2% RSD oder ≤ 0,1 SD min, je nachdem welcher Wert größer ist Bedingungen: Für das Proteinverdau-Peptid TGQAPGFSYTDANK Cytochrom C für Gradienten 30 - 90 min.
Druckbereich	Pumpendruck: 2 – 150 MPa (20 – 1500 bar, 300 – 21750 psi) 150 MPa Pumpendruck bis 1 μL/min. Bei einem Fluss über 1 μL/min ist der maximale Druck lösungsmittelabhängig und nimmt linear ab bis 70 MPa bei 100 μL/min für 80% Acetonitril in Kanal B.
Anzahl der Lösungsmittelkanäle	2
Maximales Hubvolumen	96 μL
Dwell-Volumen (Beitrag der Pumpe zum Gradientenverzögerungsvolumen des Systems)	25 nL
Entgasung	2 Kanäle für die Nadel-Waschflüssigkeiten des Autosamplers
Kommunikation	USB: 1 USB-Port (USB 2.0, Typ "B") 1 USB-Hub mit 3 Ports (USB 2.0, Typ "A") <i>E/A-Schnittstelle</i> : 2 Dig I/O-Ports (mini-DIN), jeweils mit einem digitalen Eingang und einem Relaisausgang <i>System-Interlink:</i> 2 System Interlink-Ports (RJ45-8)
Materialien im analytischen Flussweg	Keramik, DLC (diamantähnlicher Kohlenstoff), Fluorpolymere, MP35N, PEEK, Titan, UHMW-PE Informationen zur chemischen Beständigkeit der Materialien können Sie der technischen Literatur entnehmen.

Die Leistungsdaten der Pumpe sind wie folgt spezifiziert:

9.2.3 Leistungsspezifikationen des Autosamplers

Die Leistung des Autosamplers ist wie folgt spezifiziert:

Art	Spezifikation	
Injektionsmethode	Split-Loop, direkte Inje	ktion oder Trap und Elution
Anzahl der Injektionseinheiten	1	
Druckbereich	5 – 151 MPa (50 – 1517 bar, 700 – 22002 psi)	
Injektionsvolumen-Bereich	25 μL Probenschleife (Standard)	0,01 μL - 25 μL, in Schritten von minimal 0,01 μL einstellbar
	100 μL Probenschleife	0,01 μL - 100 μL , in Schritten von minimal 0,01 μL einstellbar
	(erweitert)	Hochvolumiges Trappen durch mehrere Ansaug- und Trap-Zyklen: bis zu 500 μL
Genauigkeit des Injektionsvolumens	typisch ± 0,5% mit eine	em Injektionsvolumen von 5 μ L und Wasser
Präzision des Injektionsvolumens	0,05 - 0,20 μL	<5,0%
	0,20 - 0,50 μL	<1,5%
	0,50 - 2,00 μL	<0,5%
	Über 2 μL	<0,25%
Proben-Handling mit minimalem Verlust	Aus 3 µL können mit geeigneten Probenfläschchen 2.5 µL gewonnen werden (siehe Empfehlungen zu Kappen und Verschlüssen von Probengefäßen (▶ Seite 166))	
Linearität	r > 0,9999 (Koffein in Wasser)	
Waschen der Nadel	Tauchspülen und kont unterschiedlichen Nad	inuierliches Waschen des Nadeläußeren mit 2 lel-Waschflüssigkeiten.
	Spülen der Probenschl Waschflüssigkeit aus d Probenschleife und de aus der Dosiereinheit (eife und des Nadelinneren mit starker lem Waschport (SWP). Danach Spülen der Is Nadelinneren mit schwacher Waschflüssigkeit (W).
Anzahl der Nadel-Waschflüssigkeiten	2	
Dwell-Volumen	Bei einem Probenschleifenvolumen von 25 μL:	
(Beitrag des Autosamplers zum Gradientenverzögerungsvolumen des	In Inject-Position	55 μL
Systems)	In Bypass-Position	51 nL
Verschleppung	< 0,002% mit Koffein	
Barcodeleser	Ja	
Automatisierung des Barcodelesens	Barcodelesen:	
	• Erkennung von leeren Segmenten	
	Überprüfung von Probenhalter/Mikrotiterplatte	
	Inventarverwaltung	
Probenraumthermostatisierung: Temperaturbereich	4 - 40 °C, Kühlung ≥ 23 K unterhalb der Umgebungstemperatur bei < 80% relativer Luftfeuchtigkeit	
Temperaturgenauigkeit	-2 °C und +4 °C	

Art	Spezifikation
Temperaturstabilität	±1°C
Probenkapazität	4 Segmente für Probenhalter und Mikrotiterplatten mit SBS- Flächenformat verwendbar; plus Zusatz-Probenfläschchenhalter mit einer Kapazität von (Anzahl Probenfläschchen x AD): 12 x 22,5 mm
Probenfläschchen-Kapazität	 54 x 12 mm AD Probenfläschchen 96 x 6 mm AD Probenfläschchen 96 x 7 mm AD Probenfläschchen 96 x 8 mm AD Probenfläschchen 16 x 15 mm AD Probenfläschchen 9 x 22,5 mm AD Probenfläschchen
Mikrotiterplatten-Kapazität	96- und 384-Positionen-Mikrotiterplatten mit SBS-Flächenformat
Minimal erforderliche Probe	2 μL bei 1 μL Injektionsvolumen
USB-Kommunikation	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)
Materialien im analytischen Flussweg	Titan, Keramik, PEEK, MP35N, DLC, Saphir, Fluorpolymere Hinweis: Informationen dazu, welche Hinterspülflüssigkeit zu verwenden ist, finden Sie in Füllen des Behälters für die Hinterspülflüssigkeit (> Seite 133). Informationen zur chemischen Beständigkeit der Materialien können Sie der technischen Literatur entnehmen.
Materialien im Nadelwasch-Flussweg	PP (Polypropylen), PE, FFPM (Perfluoroelastomer), FFKM, PEEK, PA (Polyamid), PK (Polyketon), TPE (thermoplastische Elastomere)
Materialien im Flussweg der Hinterspülung	Titan, Keramik, PEEK, UHMW-PE, PP, PTFE, Silikon

9.2.4 Leistungsspezifikationen des System Controllers

Die Leistungsdaten des System-Controllers sind wie folgt spezifiziert:

Art	Spezifikation
E/A-Schnittstellen:	1 LAN-Port
	2 USB-Ports (deaktiviert)
	4 System Interlink-Ports (2 sind deaktiviert)
System-Interlink	2 System Interlink-Ports (1 Paar, RJ45-8 Stecker)

9.2.5 Leistungsspezifikationen des Säulenthermostaten

Die Leistungsdaten des Säulenthermostaten sind wie folgt spezifiziert:

Art	Spezifikation
Druckbereich	Vanquish-Ventil: 50-1500 bar
	VICI-Ventil C82: 50-1034 bar
	VICI-Ventil C2, C2N: 50-345 bar
Temperaturbereich	Min.: 5 K über Raumtemperatur
	Max.: 80 °C ohne installierter Ventile
	Max.: 60 °C mit installierten Varquish-ventilen
Temperaturgenauigkeit	±0,5 °C
Temperaturstabilität	±0,05 °C
Temperaturpräzision	±0,1 °C
Aufwärmzeit	Von 35 °C auf 65 °C ±1 °C in weniger als 12 Minuten
Säulenschaltventile	Bis zu zwei Ventile:
	Vanquish-Ventil: 2 Positionen, 6 Ports, 1500 bar
	VICI-Ventil C82: 2 Positionen, 6 Ports, 1034 bar
	VICI-Ventil C2: 2 Positionen, 6 Ports, 345 bar
	VICI-Ventil C2N: 2 Positionen, 6 Ports, 345 bar
	Weitere Informationen zu erhältlichen Säulenschaltventilen finden Sie unter Optionales Zubehör () Seite 403).
USB-Kommunikation	1 USB-Port (USB 2.0, Typ "B")
	1 USB-Hub mit 3 Ports (USB 2.0, Typ "A")
E/A-Schnittstelle	2 Dig I/O-Ports (Mini-DIN), jeweils mit einem digitalen Eingang und einem Relaisausgang
System-Interlink	2 System Interlink-Ports (RJ45-8)

Art	Spezifikation
Materialien im Flussweg	Vanquish-Ventil: Keramik, DLC, Titan
	VICI-Ventil C82: Edelstahl mit inerter Beschichtung, PAEK-Komposit
	VICI-Ventil C2, C2N: PAEK, PAEK-Komposit
	Kapillaren: MP35N (weitere Informationen finden Sie bei den Vorheizer- Spezifikationen)
	Informationen zur chemischen Beständigkeit der Materialien können Sie der technischen Literatur entnehmen.

Passiver Vorheizer

Art	Spezifikation
Temperaturbereich	Wie Säulenraum
Materialien im Flussweg	MP35N
Biokompatibel	Ja
Erwärmtes Volumen	1 μL
Gesamtvolumen	6 μL
10 Zubehör, Ersatzteile und Verbrauchsmaterialien

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

10.1 Allgemeine Informationen

Das System 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

In diesem Abschnitt wird der Inhalt des Zubehörkits aufgeführt. 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 Systems finden Sie in der beiliegenden Zubehörliste.

Informationen zur Nachbestellung finden Sie unter Ersatzteile und Verbrauchsmaterialien () Seite 406).

10.2.1 System

Kapillaren

Detaillierte Informationen zur Verwendung der Kapillaren finden Sie in den Flussdiagrammen, die dem Zubehörkit beiliegen.

Artikel	Menge im Zubehör
Kapillare, nanoViper, Fused-Silica-/PEEK-ummantelt, 1/32", ID x AD x Länge 20 μm x 360 μm x 550 mm, für Nano-LC- Anwendungen	2
Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 20 μm x 280 μm x 1 m, für Nano-LC-Anwendungen	1
Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 50 μm x 360 μm x 550 mm, für Kapillar- und Mikro-LC-Anwendungen	4
Kapillare, Pumpe-Autosampler, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 50 μm x 360 μm x 350 mm, für Kapillar- und Mikro-LC-Anwendungen	1
Kapillare, Autosampler-Säule/Säulenthermostat, nanoViper, Fused-Silica-/PEEK-ummantelt, 1/32", ID x AD x Länge 50 μm x 280 μm x 1 mm, für Kapillar- und Mikro-LC-Anwendungen	1
Kapillare, Säule-Detektor, Fused Silica, ID x AD x Länge 20 μm x 280 μm x 1,5 m, für Nano-LC-Anwendungen	1
Kapillare, Säule-Detektor, Fused Silica, ID x AD x Länge 50 μm x 280 μm x 1,5 m, für Kapillar- und Mikro-LC-Anwendungen	1

Kapillarzubehör

Artikel	Menge im Zubehör
Viper-Verbindungsstück	1
Verschlussfitting, Viper, Titan	3
Rändelschraube für nanoViper-Fittinge	2

Waschsystem-Zubehör

Artikel	Menge im Zubehör
Schlauch, Silikon	3m
Schlauchverbinder, gerade (5 Stück)	1

Flüssigkeitsbehälter und Zubehör für Flüssigkeitsbehälter

Artikel	Menge im Zubehör
Einsatz für Flaschen	1
Behälter, 1 L, mit Deckel	5
Behälter, 0,25 L, mit Deckel	4
Behälterdeckel	10
Schlauchführung (Durchmesser 1,6 mm), arretiert die folgenden Schläuche in den Behälterdeckeln:	10
 Lösungsmittelschläuche (Behälter zur Pumpe) 	
 Schläuche für die Nadel-Waschflüssigkeit (Behälter zum Degaser für die Waschflüssigkeit in der Pumpe und Behälter zur Wasch- pumpe im Autosampler) 	
Schlauchführung (Durchmesser 3,0 mm), arretiert den Hinterspül- schlauch in den Behälterdeckeln	10
Verschlusskappe zum Verschließen der Behälterdeckelöffnungen (20 Stück)	2
Lösungsmittelfilter, Filterhalter (Filterfritten oder Schlauchadapter nicht enthalten)	6
Schlauchadapter	6
Lösungsmittelfilter, Filterfritte, 10 µm	10

Werkzeuge

Artikel	Menge im Zubehör
Dichtringwerkzeug	1
Spritze, Plastik, 12 mL	2
Drehmoment-Schraubendreher, Viper, Basiskomponente mit Verlängerung und Steckaufsätzen zum Festziehen der NanoViper- Verbindungen und des Autosampler-Nadelsitzes	1
Kapillarschneider für PEEK-Schläuche	1
Kapillarschneider für Fused Silica-Schläuche	1

Kabel

Artikel	Menge im Zubehör
Digitales I/O-Signalkabel, 5 m	1
LAN-Kabel, (RJ45), 5 m	1
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0, Länge: 0,5 m Zur Verbindung der Pumpe mit dem Autosampler.	1
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0, Länge: 5 m Zur Verbindung der Pumpe mit dem Computer.	1

Autosampler-Spezifische Ersatzteile

Artikel	Menge im Zubehör
Nadelsitz	1
Probenhalter, für 54 x 12 mm AD Probenfläschchen Hinweis: Die Probenhalter sind mit einem 2D-Barcode für die Vanquish-Probenhalter-Formaterkennung ausgestattet.	4
Probengefäß-Kit mit Mikrotiterplatten, Matten, Verschlussfolie, Probenfläschchen und Kappen mit vormontierten Septen	1

Sonstiges

Artikel	Menge im Zubehör
Drainage-Kit, Vanquish-System, mit: • L-Stück (4 Stück, verschiedene Größen) • Ablaufschlauch, 11,4 m Länge • Klemme (2 Stück)	1
Sicherungskit, Vanquish-System Das Kit beinhaltet die 5AT-Sicherungen für die Pumpe, den Autosampler und den Säulenthermostaten (optional).	1
Flussdiagramm	1
Cytochrom-Testprobe	1
Papieretikett für den magnetischen Etikettenrahmen	1
Spiralschlauch zum Umwickeln von Flüssigkeitsschläuchen und Kabeln	1
Halter für die Trap-Säulen-Kartusche	1
Acclaim™ PepMap™ 100 C18 HPLC-Säule (3 Stück)	1
Kappe zum Schließen von offenen Lösungsmittelschlauchenden (wenn nicht mit der Pumpe verbunden)	6
Kappe zum Schließen von offenen Ports an der Pumpe (wenn keine Lösungsmittelschläuche angeschlossen sind)	6

10.2.2 Säulenthermostat

Im Zubehörkit für den Säulenthermostat befinden sich folgende Komponenten.

Artikel	Menge im Zubehör
Säulenbefestigungsclips (6 Clips)	1
Säulen-ID-Stecker (1 Satz bestehend aus 2 Stück)	1
System-Interlink-Kabel, Länge 0,5 m Zur Verbindung des Säulenthermostaten mit dem Autosampler.	1
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0, Länge: 1 m Zur Verbindung der Pumpe mit dem Säulenthermostaten.	1

10.3 Optionales Zubehör

Probenschleifen

Beschreibung	BestNr.
Probenschleife, biokompatibel, MP35N, Volumen: 10 μ L	6252.1960
Probenschleife, biokompatibel, MP35N, Volumen: 25 μL (Standard-Probenschleife)	6252.1940
Probenschleife, biokompatibel, MP35N, Volumen: 100 μ L	6252.1950

Probenhalter und Mikrotiterplatten

Beschreibung	BestNr.
Probenhalter, für 54 x 12 mm AD Probenfläschchen	6850.1023
Nilt 2D-Barcode für die Vanquisn-Probennalter-Format-Erkennung	
Probenhalter, für 96 x 6 mm AD Probenfläschchen	6850.1026
Probenhalter, für 96 x 7 mm AD Probenfläschchen	6850.1030
Mit 2D-Barcode für die Vanquish-Probenhalter-Format-Erkennung	
Probenhalter, für 96 x 8 mm AD Probenfläschchen	6850.1034
Mit 2D-Barcode für die Vanquish-Probenhalter-Format-Erkennung	
Probenhalter, für 16 x 15 mm AD Probenfläschchen	6851.1030
Mit 2D-Barcode für die Vanquish-Probenhalter-Format-Erkennung	
Probenhalter, für 9 x 22,5 mm AD Probenfläschchen	6851.1020
Mit 2D-Barcode für die Vanquish-Probenhalter-Format-Erkennung	
Bestellinformationen zu Verbrauchsmaterialien für den Autosampler Mikrotiterplatten, Verschlussmatten, Probenfläschchen und Verschl Sie unter thermofisher.com/chromatographyconsumables.	r wie usskappen finden

Säulenthermostat-Schaltventile und -Zubehör

Artikel	BestNr.
Säulen-ID-Stecker (1 Satz bestehend aus 2 Stück)	6732.0610
Passiver Vorheizer, 0,1 mm x 680 mm, MP35N	6732.0184
Säulenschaltventil 2p-6p, niedrige Dispersion,150 MPa, biokompatibel	6250.1520
Säulenschaltventil 2p-6p, niedrige Dispersion, VICI, C82	6041.0004B
Säulenschaltventil, 2p-6p, niedrige Dispersion, 1/16", 0,25 mm, PAEK	6821.5120
Säulenschaltventil, 2p-6p, niedrige Dispersion, 1/32", 0,1 mm, PAEK	6820.6232
Verschlussfitting für Säulenschaltventil, 2p-6p, niedrige Dispersion, 1/32", 0,1 mm, PAEK (6 Stück)	6820.1320
Klemmring für VICI-Schaltventil	6253.1651

USB-Isolator

Artikel	BestNr.
USB-Isolator Der USB-Isolator verbessert die Zuverlässigkeit der USB- Verbindung zwischen dem USB-Port eines HPLC-Systemmoduls und dem Datensystemrechner.	6287.0540

System-Arbeitstische und -Stabilisatoren

Artikel	BestNr.
Bench Clamp Kit Stabilisiert Systemtürme mit einer Höhe von max. 139 cm	6036.1740
Stack Stabilizer Kit Stabilisiert Systemtürme mit einer Höhe von max. 123 cm	6036.1710
IonBench mit Montage-Kit für den Systemturm Stabilisiert Systemtürme mit einer Höhe von max. 123 cm	6036.1720

Workflows

Beschreibung	BestNr.
Tandem-Workflow-Kit für Direktinjektion, Vanquish Neo	6250.1030
Das Kit enthält alle Komponenten für den fluidischen Aufbau des Vanquish Neo Tandem-Direktinjektions-Workflows. Zusätzlich zum Kit benötigt der Workflow die folgenden Systemmodule:	
 1x Vanquish Neo-System (Standardkonfiguration) 	
 1x Vanquish Neo-Säulenthermostat mit 2x Ventilen mit niedriger Dispersion (2p-6p, 150 MPa) (BestNr. 6250.1520) 	
• 1x Vanquish Neo Binäre Pumpe	
Informationen zum Aufbau des Vanquish Neo-Tandem- Direktinjektions-Workflows finden Sie im <i>Vanquish Neo System</i> <i>User Guide</i> , Version 2.0 oder höher.	

Sonstiges

Artikel	BestNr.
Vanquish Neo Standard Application Kit	6250.1010
Kit für die Reinstallation und Requalifizierung eines bestehenden Vanquish Neo-Systems. Es enthält:	
 Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 20 μm x 280 μm, für Nano-LC- Anwendungen 	
 Kapillare, Säule-Detektor, Fused Silica, ID x AD x Länge 50 μm x 280 μm, für Kapillar- und Mikro-LC-Anwendungen 	
• Kapillar-Zubehörkit, mit:	
 Kapillare, Säule-Detektor, Fused Silica, ID x AD x Länge 20 μm x 280 μm x 1,5 m, für Nano-LC-Anwendungen 	
 Kapillare, Säule-Detektor, Fused Silica, ID x AD x Länge 50 μm x 280 μm x 1,5 m, für Kapillar- und Mikro-LC- Anwendungen 	
 Kapillare, nanoViper, Fused-Silica-/PEEK-ummantelt, 1/32", ID x AD x Länge 20 μm x 360 μm x 550 mm, für Nano-LC- Anwendungen (2 Stück) 	
 Kapillare, Pumpe-Autosampler, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 50 μm x 360 μm x 350 mm, für Kapillar- und Mikro-LC-Anwendungen 	
 Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 50 μm x 360 μm x 550 mm, für Kapillar- und Mikro-LC-Anwendungen (4 Stück) 	
 Kapillarschneider f ür Fused Silica-Schl äuche 	
 Acclaim™ PepMap™ 100 C18 HPLC-Säule (3 Stück) 	
 Halter f ür die Trap-S	
 Cytochrom-Testprobe 	

10.4 Ersatzteile und Verbrauchsmaterialien

Informationen zu Bestellnummern für Workflow-spezifische Ersatzteile finden Sie im *Vanquish Neo System User Guide*, Version 2.0 oder höher.

10.4.1 Kapillaren, Kapillarzubehör und Drainage

Kapillaren

Beschreibung	BestNr.
Kapillare, Pumpe-Autosampler, nanoViper, Fused Silica ID x Länge 20 µm x 350 mm	6250.5240
Kapillare, Autosampler-Säulenthermostat/Säule, nanoViper, Fused-Silica-/PEEK-ummantelt, 1/32", ID x Länge 20 μm x 550 mm	6250.5260
Kapillare, Pumpe-Autosampler, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x Länge 50 μm x 350 mm	6250.5540
Kapillare, Autosampler-Säulenthermostat/Säule, nanoViper, Fused-Silica-/PEEK-ummantelt, 1/32", ID x Länge 50 μm x 550 mm	6250.5560
Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 20 μm x 280 μm x 1 m, für Nano-LC-Anwendungen	6041.5292
Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 20 μm x 360 μm x 1 m, für Nano-LC-Anwendungen	6041.5293
Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 50 μm x 280 μm x 1 m, für Nano-LC-Anwendungen	6041.5294
Kapillare, Säule-Detektor, nanoViper, Fused-Silica-/PEEK- ummantelt, 1/32", ID x AD x Länge 50 μm x 360 μm x 1 m, für Nano-LC-Anwendungen	6041.5295

Kapillarzubehör

Beschreibung	BestNr.
Verschlussfitting, Viper, biokompatibel	6040.2303
Viper-Verbindungsstück	6040.2304

Kapillarwerkzeuge

Beschreibung	BestNr.
Kapillarschneider für Fused Silica-Schläuche	6720.0016
Kapillarschneider für PEEK-Schläuche	6300.0401
Drehmoment-Schraubendreher, Viper, Basiskomponente mit Adapter und Steckaufsätzen zum Festziehen der NanoViper- Verbindungen und des Autosampler-Nadelsitzes	6250.2110

Drainage-Kit

Beschreibung	BestNr.
Drainage-Kit, Vanquish-System, mit:	6036.1120
 L-Stück (4 Stück, verschiedene Größen) 	
 Ablaufschlauch, 6,25 m Länge 	
 Clip zur F ührung der Ablaufschl äuche (2 St ück) 	

10.4.2 System-Ersatzteile und Werkzeuge

System-Ersatzteile

Beschreibung	BestNr.
Solvent Rack (ohne Einsatz für Flaschen)	6036.1350
Einsatz für Flaschen im Solvent Rack	6036.1371
Vanquish Display	6036.1180
Vanquish System Controller	6036.1130-01

Die Ersatzteile für den Vanquish System Controller finden Sie unter System Controller-Ersatzteile (> Seite 408).

Werkzeuge

Beschreibung	BestNr.
Systemsockelschlüssel	6036.1160
Drehmoment-Schraubendreher, Viper, Basiskomponente mit Adapter und Steckaufsätzen zum Festziehen der NanoViper- Verbindungen und des Autosampler-Nadelsitzes	6250.2110
Werkzeugkit, VN, mit:	6250.2100
 Schraubendreher Torx T10, L-förmig 	
 Schraubendreher, Torx T20, L-förmig 	
 Schraubendreher, Torx T30 	
 Sechskantschlüssel, Größe 6 	
 Dichtringwerkzeug (Pumpe) 	
 Drehmoment-Schraubendreher, Viper, Basiskomponente mit Verlängerung und Steckaufsätzen zum Festziehen der NanoViper-Verbindungen und des Autosampler-Nadelsitzes 	
 Kapillarschneider f ür PEEK-Kapillaren 	
 Kapillarschneider f ür Fused-Silica-Kapillaren 	

10.4.3 System Controller-Ersatzteile

Beschreibung	BestNr.
SD-Karte	6439.0034
System-Interlink-Kabel, Länge: 0,5 m Zur Verbindung der Pumpe mit dem Autosampler.	6310.2012
System-Interlink-Kabel, Länge: 0,25 m Zur Verbindung folgender Komponenten: • Systemsockel zum System-Controller • System-Controller zur Pumpe	6310.2013

Netzkabel

Beschreibung	BestNr.
Netzkabel, Australien	6000.1060
Netzkabel, China	6000.1080
Netzkabel, Dänemark	6000.1070
Netzkabel, EU	6000.1000
Netzkabel, Indien, SA	6000.1090
Netzkabel, Italien	6000.1040
Netzkabel, Japan	6000.1050
Netzkabel, UK	6000.1020
Netzkabel, US	6000.1001
Netzkabel, Schweiz	6000.1030

10.4.4 Lösungsmittel und Waschsysteme

Behälter und Behälterteile

Beschreibung	BestNr.
Behälter, 1 L, mit Deckel	2270.0012
Behälter, 0,25 L, mit Deckel	2270.0026
Deckel für Behälter, Schraubdeckel (4 Stück)	6270.0013
Verschlusskappe zum Verschließen der Behälterdeckelöffnungen (20 Stück)	6000.0047
Schlauchführung (Durchmesser 1,6 mm), arretiert die folgenden Schläuche in den Behälterdeckeln (5 Stück):	6261.0216
 Lösungsmittelschläuche (Behälter zur Pumpe) 	
 Schläuche für die Nadel-Waschflüssigkeit (Behälter zum Degaser für die Waschflüssigkeit in der Pumpe und Behälter zur Wasch- pumpe im Autosampler) 	
Schlauchführung (Durchmesser 3 mm), arretiert den Hinterspül- schlauch in den Behälterdeckeln (5 Stück):	6000.0042
Verschlusskappen und Schlauchführungen (Durchmesser 1,6 mm) für Deckel der Hinterspülbehälter, Kit mit	6030.9101
 Verschlusskappe zum Verschließen der Behälterdeckelöffnungen (10 Stück) 	
 Schlauchführung für Hinterspülschlauch, arretiert den Schlauch im Behälterdeckel (5 Stück) 	

Lösungsmittelschläuche, Schläuche für die Waschflüssigkeitsbehälter und Lösungsmittelfilter

Beschreibung	BestNr.
Kit Lösungsmittelschläuche und Waschschläuche für das Waschen des Nadelinneren und Nadeläußeren, enthält die Schläuche von den Behältern und Labels für die Lösungsmittel	6250.2305
Lösungsmittelschläuche (A, B) von Lösungsmittelbehälter zum Pumpenkopf	6251.0096
Das Niederdruck-Schlauchkit für die Pumpe enthält auch die Lösungsmittelschläuche. Siehe Pumpen-Ersatzteile (▶ Seite 411).	
Lösungsmittelfilter, Filterhalter (6 Stück) (Filterfritten oder Schlauchadapter nicht enthalten)	6268.0115
Schlauchadapter (1/8-1/16) (6 Stück)	6850.1056
Lösungsmittelfilter, Filterfritte, biokompatibel, 10 μ m (10 Stück)	6268.0117

Schläuche und Teile für Waschsysteme

Beschreibung	BestNr.
Peristaltik- und Waschschlauch-Kit	6044.1150
Das Kit enthält Peristaltikschlauch (PharMed), Silikonschlauch und Schlauchverbinder für	
 Hinterspülung in Pumpe und Autosampler 	
 Drainagepumpe im Autosampler 	
ACHTUNG: Verwenden Sie den dicken Silikonschlauch (und die entsprechenden Schlauchverbinder) im Autosampler. Verwenden Sie den dünnen Silikonschlauch (und die entsprechenden Schlauchverbinder) in der Pumpe. Verwenden Sie die schraubbaren Schlauchverbinder für den Pumpenkopf.	
Detektor der Hinterspülung (Tropfendetektor) mit Trichter	6044.1898A
VN-Kit Hinterspüldichtungen und Schrauben (VN-Pumpe), Hinterspülplatte/Pumpenkopf, mit:	6251.2234
 Pumpenkopfdichtung, f ür Hintersp ülplatte (O-Ring-Dichtung, 11x1,5) (2 Dichtungen) 	
 Pumpenkopfdichtung, f ür Schraube Hintersp ülplatte (O-Ring- Dichtung, 6,5x1) (8 Dichtungen) 	
 Schraube, Hinterspülplatte (8 Schrauben) 	
Kolbendichtung (RP) und Hinterspüldichtung, Pumpenkopf (jeweils 2 Stück)	6251.1143
Nadelwaschkit, PEEK, mit	6252.1930
 Waschport-Abfallkapillare, PEEK (2 Stück) 	
 Verbindung vom Auslass der Waschpumpe zum Waschport 	

10.4.5 Pumpen-Ersatzteile

Pumpenkopf und Pumpenkopfteile

Beschreibung	BestNr.
Pumpenkopf (komplette Einheit)	6251.1102
Kolben	6251.1155
Dichtringwerkzeug	6251.5122
Kolbendichtung (RP) und Hinterspüldichtung, Pumpenkopf (jeweils 2 Stück)	6251.1143
Stützring (2 Stück)	6040.0012
Kugelventil, Einlasskugelventil	6251.1153
Kugelventil, Auslasskugelventil	6044.2310
Kapillarkit, Pumpenkopf, mit:	6251.0086
Kapillare vom Pumpenkopf zum Flowmeter	
VN-Kit Hinterspüldichtungen und Schrauben (VN-Pumpe), Hinterspülplatte/Pumpenkopf, mit:	6251.2234
 Pumpenkopfdichtung, f ür Hintersp ülplatte (O-Ring-Dichtung, 11x1,5) (2 Dichtungen) 	
 Pumpenkopfdichtung, f ür Schraube Hintersp ülplatte (O-Ring- Dichtung, 6,5x1) (8 Dichtungen) 	
 Schraube, Hinterspülplatte (8 Schrauben) 	

Flowmeter

Beschreibung	BestNr.
Filterkit, Flowmeter, mit Filterschraube, Schraubendichtung und Filterfritte (2 $\mu\text{m})$	6251.0111
Um den Filter oder die Filterfritte bei beiden Flowmeter- Eingängen zu tauschen, benötigen Sie zwei Kits.	
Flowmeter-Abfallleitung, siehe Niederdruck-Schlauchkit	

Niederdruck-Schlauchkit

Beschreibung	BestNr.
Niederdruck-Schlauchkit, Pumpe, mit: • Flowmeter-Abfallleitung vom Flowmeter zum Waste-Ablauf • Lösungsmittelschläuche (A, B) von Absperrventilen zu Pumpenköpfen	6251.0150
 Lösungsmittelschläuche vom Degaser zum Absperrventil 	
Siehe auch Lösungsmittel und Waschsysteme (> Seite 409) und Kapillaren, Kapillarzubehör und Drainage (> Seite 406).	

Wartungskit

Beschreibung	BestNr.
Wartungskit, Pumpe, mit:	6251.0130
 Schrauben, Hinterspülplatte (8 Schrauben) 	
 Silikonschlauch (durchsichtiger Schlauch) 	
 Peristaltikschlauch (PharMed, weißer Schlauch) 	
 Schlauchverbinder (gerade) 	
 Schlauchverbinder (90°-Winkel) 	
 Schlauchverbinder (ID 1/16") (4 Stück) 	
 Kolbendichtung (4 Stück) 	
 Pumpenkopfdichtung, für Hinterspülplatte (O-Ring-Dichtung, 11x1,5) (2 Stück) 	
 Pumpenkopfdichtung, f ür Schrauben Hintersp ülplatte (O-Ring- Dichtung, 6,5x1) (8 St ück) 	
 Hinterspüldichtung (4 Stück) 	
 Lösungsmittelfilter, Fritte (biokompatibel, 10 μm) (10 Stück) 	

Sonstiges

Beschreibung	BestNr.
NanoViper-Kapillare, Fused Silica, I.D. x Länge 20 μm x 350 mm, zur Verwendung als Gegendruckkapillare während der Pumpenwartung	6250.5240
Verschlussfitting, Viper, biokompatibel	6040.2303
Spritzen- und Schlauchkit, mit • Plastikspritze, 12 mL (5 Stück) • Silikonschlauch (AD x ID 2,8 x 1,3), 3 m	6000.0010
Spülschlauch, Silikon (AD x ID 2,8 x 1,3), 3 m	6007.9100
Reinigungsstäbchen (25 Stück)	6040.0007
Sicherungskit, Vanquish-System Verwenden Sie für Pumpe, Autosampler und Säulenthermostat die Sicherungen 5AT, 230 V AC, träge.	6036.0002

10.4.6 Autosampler-Ersatzteile

Kapillaren

Beschreibung	BestNr.
Verschlussfitting, Viper, biokompatibel	6040.2303
NanoViper-Kapillare, ID x L 50 µm x 120 mm	6252.5124
Abfallkapillare mit T-Stück	6252.1930
Weitere Kapillaren finden Sie unter Kapillaren, Kapillarzubehör und Drainage (▶ Seite 406).	

Flusskomponenten

Beschreibung	BestNr.
Kapillarkit für die inneren Autosampler-Flusskomponenten, bestehend aus:	6252.1920
 Viper-Kapillare, Drucksensor-Einlass 	
 Viper-Kapillare, Drucksensor-Auslass 	
 Viper-Kapillare, Dosiereinheit 	
 Viper-Kapillare, Kugelventil 	
 Waschport-Abfallkapillare, PEEK (2 Stück) 	
 NanoViper-Kapillare, Fused Silica/PEEK-ummantelt, 1/32" 	
Kopf der Dosiereinheit, 100 μL Volumen	6252.1753
Nadelsitz	6252.2470
Nadeleinheit	6252.1130
Kugelventil, einschließlich Kapillare	6252.2310
Waschsystemschläuche finden Sie in Lösungsmittel und Waschsyste	me (🕨 Seite 409).

Sonstiges

Beschreibung	BestNr.
Gerätetüren-Kit mit rechter Tür und linker Tür	6850.0100
Isolierabdeckung	6252.1647
Transportsicherung für den Drehteller	6850.5849
Drehmoment-Schraubendreher zum Festziehen der NanoViper- Verbindungen und des Nadelsitzes	6250.2110
Spülschlauch, Silikon (AD x ID 2,8 x 1,3), 3 m	6007.9100
Sicherungskit, Vanquish-System Verwenden Sie für Pumpe, Autosampler und Säulenthermostat die Sicherungen 5AT, 230 V AC, träge.	6036.0002

10.4.7 Ersatzteile des Säulenthermostaten

Beschreibung	BestNr.
Säulenbefestigungsclips (6 Clips)	6722.0290
Abdeckungen (1 Satz bestehend aus 2 Stück)	6253.1647
Gerätetüren-Kit mit rechter Tür und linker Tür	6250.1620
Rotordichtung, 2p-6p, C72/C82, E5-Ventil, 1/16", 0,15 mm	6041.0008B
Rotordichtung, 2p-6p, C2/C3-Ventil, 1/16", 0,25 mm	6820.0035
Stator, 2p-6p, C72/C82-Ventil, 1/16", 0,15 mm, N60C	6041.0007
Stator, 2p-6p, C1/C2-Ventil, 1/16", 0,25 mm	6820.0034
Sicherungskit, Vanquish-System	6036.0002
Verwenden Sie für Pumpe, Autosampler und Säulenthermostat die Sicherungen 5AT, 230 V AC, träge.	

10.4.8 Schnittstellen-Kabel und Netzkabel

Schnittstellen-Kabel

Beschreibung	BestNr.
System-Interlink-Kabel, Länge: 0,5 m Zur Verbindung der Pumpe mit dem Autosampler.	6310.2012
System-Interlink-Kabel, Länge: 0,25 m Zur Verbindung folgender Komponenten: • Systemsockel zum System-Controller • System-Controller zur Pumpe	6310.2013
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0, Länge: 0,5 m Zur Verbindung der Pumpe mit dem Autosampler.	6720.8910A
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0, Länge: 1 m Zur Verbindung der Pumpe mit dem Säulenthermostaten	6035.9035A
USB-Kabel, Typ A auf Typ B, High-speed, USB 2.0, Länge: 5 m Zur Verbindung der Pumpe mit dem Computer.	6911.0002A
Digitales I/O-Signalkabel, 6-polig, Kabellänge: 5 m	6036.0006

Netzkabel

Beschreibung	BestNr.
Netzkabel, Australien	6000.1060
Netzkabel, China	6000.1080
Netzkabel, Dänemark	6000.1070
Netzkabel, EU	6000.1000
Netzkabel, Indien, SA	6000.1090
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 und zur Verwendung der digitalen E/A-Schnittstellen.

11.1 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.

11.1.1 Konformitätserklärungen

CE-Konformitätserklärung

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

EAC-Konformitätserklärung

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

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 Richtlinien 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

UKCA-Konformitätserklärung

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

Einhaltung der UL/CSA 61010-1-Richtlinien

Das NRTL-Zeichen auf dem Gerät (zum Beispiel cTUVus-Zeichen oder CSA-Zeichen) zeigt an, dass das Gerät die Anforderungen geltender Normen erfüllt.

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 232: 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	VN-A10-A-02, VN-C10-A-01, VN-P10-A-01, VN-S10-A-01
1.0	VN-A10-A-02, VN-C10-A-01, VN-P10-A-01, VN-S10-A-01

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

11.2 Digital I/O (Anschlussbelegung)

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

- einen digitalen Eingang
- einen Relaisausgang



Abbildung 233: Digital I/O-Port

Pin	Beschreibung - Signalname
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	 Digitaleingang; das Referenzpotential ist Ground. Beachten Sie folgende Hinweise: Die maximale Eingangsspannung darf am Eingang +5 V in Bezug auf Ground nicht überschreiten. Die minimale Eingangsspannung darf das Groundpotential nicht unterschreiten.
5	Weiß	Relaisausgang — Relay_COM	Potentialfrei	Mittelkontakt für NO und NC
6	Braun	Relaisausgang — Relay_NO	Potentialfrei 0-24 V, 0-100 mA	Schließer

Sehen Sie dazu auch

Anschließen eines Massenspektrometers (> Seite 117)

Index

Α

Abfallbehälter 125
Abmessungen
Additive
allgemeine Kompatibilität 144
Verwenden 145
Algen 145
Ändern
Hinterspülflüssigkeit 170
Nadel-Waschflüssigkeit 170
Anforderungen
Abstände vom System 79
Arbeitstisch
Netz
Netzkabel 80
Stromversorgung (Hinweise)
Anforderungen an den Arbeitstisch 78
Anforderungen an den Aufstellungsort 78
Betriebsbedingungen 81
Kondensation 81, 115, 314
Stromversorgung (Hinweise)
Anschließen der Netzkabel 115
Anschluss
Netzkabel 314
Anschlüsse
System Controller 108, 312
Systemmodule 107
Anschlüsse der Systemmodule 107
Äquilibrierdauer 174
Auslasskugelventil
Reinigen 227
Tauschen 226
Außerbetriebnahme
kurzzeitig 175
langfristig 175
Autosampler
Besonderheiten 35
Funktionsprinzip 46
Innenansicht 44
Inventory Scan (Bestandsprüfung) 148
Moduleinschub entfernen 331
Thermostatisierung 173
Transportsicherungen
Vorbereiten für das Entfernen
Autosampler-Ventile

Port-Belegung	. 48
Undichtigkeit	296

В

Barcode (Vanquish)	51, 165
Behälter	
Kolbenhinterspülung	133
Lösungsmittel	128
Waschen der Nadel	130
Benachrichtigungsfenster	344
Betrieb	141
allgemeine Hinweise	148
Bedienelemente	150
Chromatographie-Datensystem	64
Sicherheitshinweise	142
Unterbrechung (kurz)	175
Unterbrechung (lang)	175
Vanquish User Interface	64
Betriebsbedingungen	81
Belüftung	81
Elektromagnetische Interferenz	82
Höhe	388
Luftfeuchtigkeit	81, 388
mechanische Stöße	82
Temperatur	81, 388
Vibrationen	82

С

CE-Kennzeichnung	418
Chloridkonzentration	145
cTUVus-Zeichen	419

D

Dekontaminierung	190
Detaillierter Dichtigkeitstest (Pumpe)	386
Dichtigkeitstest	
Drainagepumpen-Schlauch	265
Hinterspülung	279
Systemdichtigkeit	385
Dichtungsband	166
Dig I/O	421
Digital I/O 107,	421

Display

Abbauen	317
Anpassen	158
Anschließen	111
Installieren	100
Positionen	158
seitliche Position einstellen	159
Tauschen	317
Wartung	309
Displayposition anpassen	158
Drainage	123
Drainagepumpen-Schlauch	265
Tauschen	267
Test auf Dichtigkeit	266
Testen	265
Testen auf Blockierungen	266
Drehmoment-Schraubendreher	121
Drehteller	49, 50
Beladen	163
Farbkennzeichnung	50
rotieren	164
Zusatz-Fläschchenhalter	50
Zusatz-Flaschchenhalter	50

Ε

EAC-Kennzeichnung 418
Ein- und Ausschalten 62
Einhaltung der UL/CSA-Richtlinien 419
Einlasskugelventil
Reinigen 227
Tauschen 224
Einsatz für Flaschen
Einsetzen 323
Entfernen 322
Einschalten des Systems 149
Einschalten des Systems nach der Installation
Einzuhaltende Abstände vom System
Emissionsschalldruckpegel 389
Ersatzteile 406

F

FCC	419
Fehlersuche	343
Allgemeine Informationen	344
Basislinienprobleme	379
Druckprobleme	376
instabile Analyse	375

Längere Ladezeit	380
Leak-Alarm	385
Meldungen	346
Negative Peaks	373
Peakform-Probleme	372
Peakprobleme 371,	373
Probleme bei der Stromversorgung	381
Probleme mit der Kommunikation	382
Retentionszeitschwankungen	377
Schlechte Präzision der Peakflächen	374
schlechte Reproduzierbarkeit	375
Skripte	369
Spikes	374
Störpeaks	373
Temperaturregelungsprobleme beim	
Autosampler	381
Temperaturregelungsprobleme beim	
Säulenthermostaten	380
Flasche	
Kolbenhinterspülung	133
Lösungsmittel	128
Waschen der Nadel	130
Flowmeter	
Abfallleitung tauschen	231
Filter tauschen	229
Wartung	229
Flussrate (Abhängigkeit mit Pumpendruck)	167
Flusssensor (Nullpunktabgleich durchführen)	171
Flussverbindungen	
Drainagepumpen-Schlauch	265
Systemdrainage	123
Waste-Leitungen	123
Führungsloch	119
Funktion "Catch Valve" 91,	305
Funktionsprinzip	
Autosampler	. 46
Pumpe	. 42
Säulenthermostat	. 54

G

Geräte-LAN
Integrieren des Systems 137
Gerätetür
Aushängen 325
Öffnen 85
Öffnungsmechanismus des Autosamplers 85
Schließen 85
Schließmechanismus des Autosamplers 85
Tauschen 325
Gesetzliche Bestimmungen 32
Gewicht 389
Griffe
Befestigen am System 340

Н

Handschuhe	28
Hauptnetzschalter	62, 135
Hinterspüldetektor (Tausch)	293
Hinterspüldichtung	
Entfernen	215
Installieren	217
Tauschen	215
Hinterspülflüssigkeit	
Ändern	170
Hinterspülung	60, 279
Auffüllen	133
Hinterspüldetektor	293
Hinterspül-Einlassschlauch	284
Hinterspülflüssigkeit	133
Hinterspülpumpe	289
Hinterspülpumpen-Einlassschlauch	287
Hinterspülschlauch zum Behälter	281
Hinterspülschläuche in der Pumpe	292
Peristaltikschlauch	289
Spülen	136
Test auf Dichtigkeit	279
Wartung	279
Hinweise	
Betrieb	142
Installation	76
Wartung	179
-	

I

Inject-Response	117
Injektionsventil	

Port-Belegung 48
Undichtigkeit 296
Innenansicht
Autosampler 44
Pumpe 41
Säulenthermostat 53
Installation
Anforderungen an den Aufstellungsort 78
Anschließen der Netzkabel 115
Kapillaren und Schläuche 119
Sicherheitshinweise
System 83
Waste-Leitungen 123
Interlink 39, 107
Inventory Scan (Bestandsprüfung) 148
Isolierabdeckung
Entfernen 98
Installieren 99

J

Jährliche War	rtungsarbeiten	187
---------------	----------------	-----

Κ

Kabelhalterung	116
Kalibrieren kundenspezifischer Lösungsmittel	
	168
Kapillaren	
Führung	119
Installieren	120
Kolben	
Entfernen	200
Installieren	201
Reinigen	202
Tauschen	200
Kolbendichtung	
Einsetzen	211
Entfernen	208
Tauschen	208
Kolbendichtungshinterspülung	. 60
Kolbenhinterspülung	. 60
Kondensation	314
Konnektivitätseinstellungen	157
Kopf der Dosiereinheit	
Anschließen von Kapillaren	261
Entfernen von Kapillaren	261
Tauschen	261

Kugelventil	
Pumpenkopf	224
Waschen der Nadel	278
Kugelventil (Pumpenkopf)	
Auslassventil tauschen	226
Einlasskugelventil tauschen	224
Reinigen	227
Kurzzeitige Außerbetriebnahme	175

L

Lagerungsbedingungen	
Luftfeuchtigkeit	388
Temperatur	388
Langfristige Außerbetriebnahme	175
LAN-Kabel	
Anschließen	115
Leakerkennung	. 57
Leckage-Flüssigkeit	123
LED-Leiste 150, 153,	344
Leistung optimieren	167
Lieferumfang	. 69
Lösungsmittel	
allgemeine Kompatibilität	144
Chloridkonzentration	145
Kalibrierung	168
pH-Bereich	144
Pufferkonzentration	145
Qualität und Filtration	145
Verwenden	145
Wechseln	146
Lösungsmittelbehälter 128,	146
Lösungsmittelfilter	
Filter tauschen	240
Fritte tauschen	240
Wartung	234
Zusammensetzen 128,	241
Lösungsmittelschläuche	
Absperrventil zum Pumpenkopf	239
Behälter zu Pumpe	234
entleeren	234
Wartung	234
Lösungsmitteltyp	
Ändern 162,	170
Lösungsmitteltyp ändern 162,	170
Lösungsmittelventil	
Port-Belegung	. 49
Undichtigkeit	296

Μ

Massenspektrometer (anschließen)	117
Meldungen	
allgemein	346
Autosampler	363
Benachrichtigungsfenster	344
Pumpe	352
Säulenthermostat	368
Übergreifende Meldungen	346
Mikrotiterplatte	. 50
Abdeckungen	166
Ausrichtung	164
Barcode	. 51
Moduleinschub	
Installieren	336
Zurückschicken	335
Moduleinschub (Pumpe)	
Entfernen	328

Ν

Nadeleinheit	
entsperren	249
Service-Position	248
Sichern	248
Nadel-Waschflüssigkeit	130
Ändern	170
Nadelwasch-Kugelventil	
Tauschen	278
Nadel-Waschsystem	
Auffüllen	130
Purgen	132
Schläuche tauschen 270, 272,	273
Schläuche vom Behälter 270,	273
Waschen der Nadel	172
Waschpumpen-Auslassschlauch	272
nanoViper-Kapillaren	
Anschließen	121
Netz	
Anforderungen	388
Einschalten 135,	149
Schalter 62,	135
Netzkabel 80,	314

Ρ

Passiver Vorheizer
Installieren 92
Tauschen 307
Peristaltikschlauch
Tauschen 290
pH-Bereich144
Probenhalter 50
Ausrichtung164
Barcode
Probenhalterformat-Erkennung 51
Probenhalter-Format-Erkennung 165
Probenraum 49
Drehteller 50
Thermostatisierung 173
Probenraum-Thermostatisierung 50
Puffer
Konzentration145
Verwenden 145
Pumpe
Besonderheiten 34
Funktionsprinzip 42
Innenansicht 41
Moduleinschub entfernen 328
Vorbereiten für das Entfernen
Pumpe (spülen) 187, 189
Pumpendruck (Abhängigkeit mit Flussrate) 167
Pumpenkopf
Auslasskugelventil 226
Einlasskugelventil 224
Einsetzen 197
Entfernen 195
Kugelventil 224, 227
Tauschen 195
Übersicht Teile 194
Wartung 193
Pumpenkopfdichtung
Entfernen 219
Installieren 221
Tauschen 219
Pumpen-Meldungen 352
Purgen 65

R

Ready Temp Delta	174
Ready-Zustand	174
Regelmäßige Wartungsarbeiten	186

Reinigen	190
Reparatur	183
RoHS-Kennzeichnung	418
ROTATE (Taste)	164
Rotordichtung des VICI-Schaltventils	
Tauschen	305

S

Säulen	
Installieren	94
Tauschen	298
Säulenschaltventil	54
Undichtigkeit	299
Säulenthermostat	
Besonderheiten	35
Funktionsprinzip	54
Innenansicht	53
Installieren	86
Moduleinschub entfernen	333
Thermostatisierung	173
Vorbereiten für das Entfernen	332
Säulenthermostat auspacken	72
Schläuche	120
Schlauchführungen	119
Schlauchkanal	119
Schlauchklammer	119
Schutzbrille	28
Schutzklasse	26
Schutzkleidung	28
schwaches Waschen	130
SD-Karte	
Tauschen	315
Septen	166
Seriennummernaufkleber des Systems	23
Service	177
Sicherheitshinweise	
allgemein	26
allgemeine Risiken	29
Betrieb	142
elektrische Sicherheit	28
Installation	76
Netzkabel	80
Qualifikation des Personals	27
Schutzausrüstung	27
Service	179
Verhalten im Notfall	31
Wartung	179
Sicherheitssymbole 2	2, 23

Sicherungen	324
Signalworter	22
Skript Auf System-Werkseinstellungen	270
zurucksetzen	
Skript Autostart	138, 160
Skript System reinigen	
Skript Systemdichtigkeit testen	369
Skript Systemeinrichtung starten	370
Skript Systemrückdruck testen	369
Software-Lizenzen	157
Solvent Rack	
Einsatz für Flaschen einsetzen	323
Einsatz für Flaschen entfernen	322
Tauschen	319
Spezifikationen	387
Abmessungen	389
Autosamplerleistung	392
Emissionsschalldruckpegel	389
Gewicht	389
Leistung des System-Controllers	394
Leistungsaufnahme	388
Leistungsdaten des Säulenthermosta	ten 394
Pumpenleistung	391
Systemleistung	390
Split-Loop-Prinzip	46
Spracheinstellungen	157
Spülen	
nach Betrieb	146
Pumpe	187, 189
starkes Waschen	130
Stator des VICI-Schaltventils	
Tauschen	305
Statusanzeige	
LED-Leiste	150, 344
Status-LED	150, 344
Status-LED	150, 344
Stromversorgung (Hinweise)	79, 80
Stützring	
Entfernen	203
Installieren	205
Tauschen	203
Symbole auf dem Vanquish User Interfa	ce
· · · · · · · · · · · · · · · · · · ·	155, 156
System	
äquilibrieren	162
Außerbetriebnahme	175
Beschreibung	37
Besonderheiten	

Betrieb vorbereiten	160
Erstinbetriebnahme vorbereiten 13	8, 160
Fehlersuche	369
Leistung optimieren	167
Status	155
Übersicht	37
Vorbereiten für den Transport	338
System auspacken	70
System Controller	40
Anschließen 10	8, 312
Datum	155
Entfernen	309
Firmware-Version	155
Installieren	310
Seriennummer	155
Uhrzeit	155
Wartung	309
System Controller-Anschlüsse 10	8, 312
System einschalten	
nach der Installation	135
regelmäßig	149
System-Controller	
Tausch der SD-Karte	315
System-Fehlersuche	369
System-Interlink	9, 107
System-Interlink-Kabel	
Anschließen	113
Systemsockel	38
entsperren	73
Sperren	118
Systemsockel entsperren	73
Systemsockelschloss	39

Т

150 151
151
152
151
389
388
388
389
389
388
388
388

Temperatur (Lagerung)	388
Tragegriffe	
Befestigen am System	340
Transport	
Vorbereiten des Systems	338
Transportsicherungsschraube	
Festziehen	248
Lösen	97, 249
Tropfendetektor (Tausch)	293
Tür	
Aushängen	178
Entfernen	178
Typenschild	24
Typenschilder	23

U

Überblick (Funktionen)	33
Übersicht (Systemmodule)	37
UKCA-Kennzeichnung	. 419
Undichtigkeit	. 385
Undichtigkeiten	. 385
Universal Serial Bus	. 107
USB	. 107
USB-Kabel	
Anschließen	. 111

V

Vanquish User Interface	
DHCP-Option	137
Konnektivitätseinstellungen	157
Konnektivitätsoption	137
Option Benutzerdefiniert	137
Skripte	369
Software-Lizenzen	157
Spracheinstellungen	157
Symbole	155, 156
Zeitzone	157
Vanquish-Schaltventil	55
Installieren	88
Tauschen	300
Versand	
Moduleinschub	335
Verschlüsse	166
Verschlussmatten	166
Verwendungszweck	25
VICI-Schaltventil	55
Installieren	90

W

Wartung 177,	184
Allgemeine Regeln	182
Dekontaminierung	190
Display	309
Einführung	178
Reinigen	190
Sicherheitshinweise	179
Sicherungen	324
System Controller	309
Test auf Dichtigkeit	385
Türen	325
Wartungszeitplan	184
Jährlich	187
regelmäßig	186
Wöchentlich	185
Waste-Leitungen	123
WEEE	419
werksseitige Standardwerte	
Reset	370
werksseitige Standardwerte zurücksetzen	370
Wöchentliche Wartungsarbeiten	185

Ζ

Zeitzone	157
Zubehörkit	399
Zurückschicken	
Moduleinschub	335
System	183

www.thermofisher.com

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

