



# Thermo Scientific Dionex UltiMate 3000-Serie

# WPS-3000 Autosampler (SL, TBSL, RS, TBRS, TXRS)

Bedienungsanleitung (Originalbedienungsanleitung)



Version: 1.8 Datum: September 2013

© 2013 Thermo Fisher Scientific Dok.-Nr. 4828.2201

# CE

# EG KONFORMITÄTSERKLÄRUNG

(Original-Konformitätserklärung)

Geräteart:	Thermo Scientific Dionex UltiMate 3000 - Autosampler

Typen:WPS-3000SL und WPS-3000T(B)SLWPS-3000RS, WPS-3000T(B)RS und WPS-3000TXRS

Die Dionex Softron GmbH bescheinigt hiermit, dass die oben beschriebenen Produkte den entsprechenden Anforderungen der folgenden Richtlinien entsprechen:

- Niederspannungsrichtlinie 2006/95/EG
- EMV Richtlinie 2004/108/EG

Zur Beurteilung der Erzeugnisse hinsichtlich der elektrischen Sicherheit wurde folgende Norm herangezogen:

• DIN EN 61010-1:2010 Sicherheitsbestimmungen für elektrische Mess-, Steuer-, Regel- und Laborgeräte, Teil 1: Allgemeine Anforderungen

Zur Beurteilung der Erzeugnisse hinsichtlich der Elektromagnetischen Verträglichkeit (EMV) wurde folgende Norm herangezogen:

• DIN EN 61326:2006 Elektrische Betriebsmittel für Leittechnik und Laboreinsatz EMV-Anforderungen

Diese Erklärung wird verantwortlich für den Hersteller

Dionex Softron GmbH Part of Thermo Fisher Scientific Inc. Dornierstraße 4 D-82110 Germering

abgegeben durch den Managing Director, Rüdiger Obst und den Vice President HPLC, Fraser McLeod.

Germering, den 02.09.2013

# Inhaltsverzeichnis

1	Ei	inführung	1
	1.1	Über die Bedienungsanleitung	1
	1.2	Sicherheitsinformationen	3
	1.2	2.1 Symbole am Autosampler und in der Bedienungsanleitung	3
	1.2	2.2 Sicherheitsmaßnahmen	4
	1.3	Verwendungszweck des Autosamplers	8
2	Ül	berblick	11
	2.1	Kurzbeschreibung	11
	2.2	Funktionsprinzip	12
	2.3	Autosamplerkonfigurationen	14
	2.3	3.1 Übersicht	14
	2.3	3.2 Autosampler mit Probenthermostatisierung	16
	2.3	3.3 Kombinationsmöglichkeiten von Spritzen und Schleifen	17
	2.4	Innenansicht	18
	2.5	Gerätevorderseite	20
	2.6	Geräterückseite	21
	2.6	6.1 Netzschalter	22
	2.6	6.2 Sicherungsschlitten	22
	2.6	6.3 USB-Anschluss (USB 1.1)	22
	2.6	6.4 Digital I/O	22
	2.7	Drehteller, Probenhalter und Mikrotiterplatten	23
	2.8	Probenraumbeleuchtung	25
	2.9	Fluidische Anschlüsse	26
	2.10	Leaksensor	26
	2.11	Injektionsventil	27
	2.12	Steuerung über Chromeleon	28
	2.13	Wellness, Predictive Performance und Diagnose	29
3	In	stallation	31
	3.1	Anforderungen an den Standort	31
	3.2	Auspacken	32
	3.3	Position des Autosamplers im UltiMate 3000-System	33
	3.4	Verbinden des Autosamplers	35
	3.4	4.1 Allgemeine Informationen	35
	3.4	4.2 Anschluss des USB-Kabels	35
	3.4	4.3 Anschluss des Netzkabels	36
	3.4	4.4 Anschluss des Signalkabels (Digital I/O)	36
	3.5	Einrichten des Autosamplers in Chromeleon	37
	3.5	5.1 Laden des USB-Treibers	37
	3.5	5.2 Installieren des Autosamplers	39
	3.5	5.3 Konfigurieren des Autosamplers	40
	3.6	Einrichten des Autosamplers in DCMSLink	46

4	Vorl	oereitung für den Betrieb (Inbetriebnahme)	47
	4.1 Ü	bersicht	47
	4.2 H	inweise zum Anschluss von Kapillaren	48
	4.3 A	nschließen der Waschflüssigkeit	50
	4.3.1	Anschluss der Waschflüssigkeit bei Online-Entgasung	50
	4.3.2	Anschluss der Waschflasche	52
	4.4 A	nschlüsse am Spritzenventil	53
	4.5 A	nschließen der Drainage	
	4.6 E	ntfernen von Luftblasen aus der Spritze (Prime)	
	4.7 E	insetzen und Bestücken eines Probenhalters	
	4.8 B	estücken der 5-Positionen Halter	
	4.9 A	quilibrieren des Systems	60
5	Betr	ieb und Wartung	63
	5.1 E	inschalten des Autosamplers	63
	5.2 S	atusanzeige	64
	5.3 P	robenposition	65
	5.4 S	teuerung über Chromeleon	67
	5.4.1	Verbinden mit Chromeleon	67
	5.4.2	Direkte Steuerung	68
	5.4.3	Automatische Steuerung	71
	5.5 F	unktionstasten und Menüs am Autosamplerdisplay	73
	5.5.1	Einblenden der Funktionstasten	
	5.5.2	Autosampler-Menus	
	3.6 E	Instellungen für den Autosampler-Betrieb	81
	5.0.1	Symphronigiaran von Inightionshafahl und Kalbanhüban	81 82
	5.6.2	Finschalten der Probenthermostatisierung	
	5.0.5	Ändern der Drehtellerkonfiguration	83 84
	565	Festlegen der Injektionsnarameter und Starten der Injektion	
	5.6.6	Frkennen von Undichtigkeiten (Leakerkennung)	 86
	567	Annassen von Helligkeit und Kontrast der Displayanzeige	86
	568	SmartStartup und SmartShutdown	87
	5.6.9	Empfehlungen für den Betrieb	
	5.7 S	pezielle Funktionen in Chromeleon	
	5.7.1	Aktive Überwachung von Verschleißteilen (Predictive Performance)	
	5.7.2	Diagnosetests	90
	5.7.3	Definieren der Nadelhöhe (Sample Height, Reagent Liquid Height)	91
	5.7.4	Bewegen des Drehtellers vor dem Aufziehen der Probe (Tray Shake)	94
	5.7.5	Belüften des Probenfläschchens	94
	5.7.6	Operational Qualification und Performance Qualification	94
	5.7.7	Automatischer Tray-Test	94
	5.7.8	Automatische Probengefäßerkennung	95
	5.7.9	Benutzerdefinierte Programme	96
	5.8 A	ußerbetriebnehmen des Autosamplers	98
	5.9 W	Vartung und Wartungsintervalle	100

6	Fehlersuche	103
	6.1 Übersicht	103
	6.2 Meldungen auf dem Autosamplerdisplay	104
	6.3 Diagnose-Meldungen in Chromeleon	110
	6.4 Mögliche Störungen	112
7	Service	119
	7.1 Allgemeine Hinweise und Sicherheitsmaßnahmen	119
	7.2 Beseitigen von Undichtigkeiten	120
	7.3 Wechseln der Spritze	121
	7.4 Wechseln der Pufferschleife	124
	7.5 Probennadel	125
	7.5.1 Wechseln der Probennadel	125
	7.5.2 Waschen der Probennadel	130
	7.6 Wechseln der Probenschleife	131
	7.6.1 WPS-3000SL Analytical, WPS-3000RS/TBRS/TXRS	131
	7.6.2 WPS-3000TBSL Analytical	134
	7.6.3 WPS-3000SL Semiprep	136
	7.7 Nadelsitz	139
	7.7.1 Wechseln des Nadelsitzes	139
	7.7.2 Wechseln der Nadelsitzkapillare	142
	7.8 Injektionsventil	148
	7.8.1 Prüfen des Injektionsventils auf Undichtigkeit	149
	7.8.2 Wechseln des Ventil-Pods	150
	7.8.3 Wechseln des Ventilstators	154
	7.8.4 Rotor Seal und Stator Face Seal	155
	7.9 Wechseln der Sicherungen	158
	7.10 Aktualisieren der Autosampler-Firmware	159
8	Optimieren für spezielle Anwendungen	161
9	Technische Daten	163
	9.1 WPS-3000SL Analytical und WPS-3000SL Semiprep	163
	9.2 WPS-3000TBSL Analytical	165
	9.3 WPS-3000RS	167
	9.4 WPS-3000TBRS	169
	9.5 WPS-3000TXRS	171
1	Zubehör, Ersatzteile und Verbrauchsmaterialien	173
	10.1 Standardzubehör	173
	10.2 Optionales Zubehör	178
	10.3 Ersatzteile und Verbrauchsmaterialien	183
1	Anhang - Anschlussbelegung Digital I/O	191
12	2 Index	193

# 1 Einführung

# 1.1 Über die Bedienungsanleitung

Dieses Handbuch soll Ihnen den gezielten Zugriff auf diejenigen Abschnitte ermöglichen, die Sie für den Gebrauch Ihres Thermo Scientific<sup>TM</sup> Dionex<sup>TM</sup> WPS-3000 Autosamplers benötigen. Dennoch sollten Sie, bevor Sie mit dem Autosampler arbeiten, die gesamte Anleitung einmal gründlich durchlesen, um sich einen Überblick zu verschaffen.

Die Beschreibungen innerhalb des Handbuchs beziehen sich auf die folgenden Modelle der UltiMate™ 3000-Autosamplerserie:

- WPS-3000U(T)SL Analytical und WPS-3000(T)SL Semiprep
- WPS-3000TBSL Analytical
- WPS-3000(T)RS
- WPS-3000TBRS
- WPS-3000TXRS

Für die Beschreibungen innerhalb dieses Handbuches gelten die folgenden Konventionen:

- Für die Beschreibungen wird der Ausdruck "das Gerät" oder "der Autosampler" verwendet. Bezieht sich eine Beschreibung nur auf ein bestimmtes Modell oder eine bestimmte Version ist dies entsprechend gekennzeichnet.
- Wenn nicht anders angegeben, gelten alle Beschreibungen für
  - WPS-3000SL Analytical, WPS-3000SL Semiprep und WPS-3000RS auch f
    ür die Versionen mit Probenthermostatisierung (WPS-3000TSL Analytical, WPS 3000TSL Semiprep, WPS-3000TRS).
  - ◆ Viper<sup>™</sup>-Kapillarverbindungen ebenso für nanoViper<sup>™</sup>- und gegebenenfalls andere Viper-Kapillarverbindungen.
- Die Geräteausstattung kann je nach Geräteversion (z.B. Temperaturregelung für Probentemperierung) variieren. Daher müssen nicht alle Beschreibungen auf das ausgelieferte Gerät zutreffen.
- Die optische Ausführung einzelner Bauteile kann gegebenenfalls von den Abbildungen im Handbuch abweichen. Dies hat jedoch keinen Einfluss auf die Beschreibungen.
- Die Beschreibungen in dieser Anleitung beziehen sich auf die Firmware-Version 4.13 und Chromeleon<sup>™</sup> 6.80 Service Release 13. Wenn Sie den Autosampler unter Chromeleon 7 betreiben möchten, beachten Sie die Hinweise auf Seite 28.

Das vorliegende Handbuch wurde nach bestem Wissen und Gewissen erstellt. Alle technischen Angaben und Programme wurden mit größter Sorgfalt erarbeitet. Trotzdem sind Fehler nicht ganz auszuschließen. Wir möchten deshalb darauf hinweisen, dass weder eine Garantie noch irgendeine Haftung für Folgen, die auf fehlerhafte Angaben zurückgehen, übernommen werden kann. Hinweise auf eventuelle Fehler sind jederzeit willkommen.

Die in diesem Dokument enthaltenen Angaben und Daten können jederzeit ohne vorherige Ankündigung geändert werden.

Alle Rechte vorbehalten, auch die der fotomechanischen Wiedergabe und der Speicherung auf elektronischen Medien. Kein Teil dieser Unterlagen darf in irgendeiner Form (durch Fotokopie, Mikrofilm oder andere Verfahren) ohne die schriftliche Genehmigung seitens Thermo Fisher Scientific für irgendeinen Zweck reproduziert oder unter Verwendung elektronischer Systeme verarbeitet, vervielfältigt, übertragen oder auf andere Art und Weise verbreitet werden. Dies ist unabhängig davon, auf welche Art und Weise oder mit welchen Mitteln, elektronisch oder mechanisch, dies geschieht.

#### Warenzeichen

Analyst ist ein eingetragenes Warenzeichen von AB Sciex.

Compass und Hystar sind Warenzeichen von Bruker Daltonics.

Empower ist ein Warenzeichen von Waters Corp.

MP35N ist ein eingetragenes Warenzeichen von SPS Technologies.

PEEK ist ein Warenzeichen von Victrex PLC.

Vespel ist ein eingetragenes Warenzeichen von E.I. du Pont de Nemours & Company.

Microsoft, Windows und Windows Vista sind eingetragene Warenzeichen von Microsoft Corp.

Alle anderen Warenzeichen sind Eigentum von Thermo Fisher Scientific Inc. und ihren Tochtergesellschaften.

# 1.2 Sicherheitsinformationen

Die CE- und cTUVus-Zeichen auf der Geräterückseite geben an, dass das Gerät die entsprechenden Standards erfüllt.

#### 1.2.1 Symbole am Autosampler und in der Bedienungsanleitung

Die nachfolgende Tabelle gibt einen Überblick über die auf dem Gerät verwendeten Symbole:

Symbol	Beschreibung
~	Wechselstrom
- 0	Stromversorgung eingeschaltet (–) bzw. ausgeschaltet ( <b>O</b> )
	Handverletzungsgefahr
<b>2 ()</b>	Kennzeichnung entsprechend der Richtlinie "Measures for Administration of the Pollution Control of Electronic Information Products" (China-RoHS)
	Lesen Sie im Handbuch nach, um ein Verletzungsrisiko auszuschließen bzw. Schäden am Gerät zu vermeiden.
	WEEE-Kennzeichnung (Waste Electrical and Electronic Equipment) - Weitere Informationen finden Sie unter WEEE-Information im Ordner "Installation and Qualification Documents for Chromatography Instruments".

Innerhalb des Handbuchs machen folgende Symbole auf besonders wichtige Informationen aufmerksam:

i	Hinweis:	Hier finden Sie allgemeine Informationen sowie Informationen, die Ihnen zu optimalen Ergebnissen verhelfen sollen.
⚠	Vorsicht:	Falls Sie diese Informationen ignorieren, kann dies zu falschen Ergebnissen oder zu Schäden am Gerät führen.
STOP	Warnung:	Wenn Sie diese Informationen ignorieren, schaden Sie möglicherweise Ihrer Gesundheit.

#### 1.2.2 Sicherheitsmaßnahmen

Wenn Sie mit analytischen Geräten arbeiten, müssen Sie die Gefahren kennen, die beim Umgang mit chemischen Stoffen auftreten können.

**Hinweis:** Bevor Sie mit dem Autosampler zu arbeiten beginnen, lesen Sie diese Anleitung vollständig, so dass Sie mit dem Inhalt vertraut sind.

Warnung: Alle Benutzer des Gerätes müssen die folgenden Sicherheitshinweise und alle weiteren Sicherheitshinweise in dieser Anleitung beachten, um bei Betrieb, Wartung und Service eine Gefährdung ihrer Person oder Schäden am Gerät zu vermeiden.

> Beachten Sie etwaige Warnaufkleber auf dem Gerät und die Informationen in den entsprechenden Kapiteln der *Bedienungsanleitung*.

#### Schutzausrüstung

Tragen Sie bei allen Arbeiten an und in der Nähe des HPLC-Systems persönliche Schutzausrüstung (Schutzkleidung, Sicherheitshandschuhe, Schutzbrille), die der Gefährdung durch die mobilen Phase und Probe entspricht. Informationen zum richtigen Umgang mit konkreten Substanzen und Empfehlungen für konkrete Gefahrensituationen entnehmen Sie bitte dem Sicherheitsdatenblatt der Substanzen, mit denen Sie umgehen. Beachten Sie die Richtlinien der Guten Laborpraxis (GLP).

In der Nähe Ihres Arbeitsplatzes sollten sich auch 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.

#### • Gefährliche Substanzen

Viele organische Lösungsmittel, mobile Phasen und Proben sind gesundheitsschädlich. Vergewissern Sie sich, dass Sie die toxischen und infektiösen Eigenschaften der von Ihnen eingesetzten Substanzen kennen. Bei vielen Substanzen sind Ihnen deren toxische oder infektiöse Eigenschaften eventuell nicht bekannt. Behandeln Sie Substanzen im Zweifelsfall, als würden sie eine gesundheitsschädliche Substanz enthalten. Anweisungen zum richtigen Umgang mit konkreten Substanzen entnehmen Sie bitte dem Sicherheitsdatenblatt (SDB) des jeweiligen Herstellers. Beachten Sie die Richtlinien der Guten Laborpraxis (GLP).

Entsorgen Sie Abfälle der Substanzen umweltgerecht und entsprechend den lokalen Bestimmungen. Vermeiden Sie die Ansammlung von entzündlichen, toxischen und/oder infektiösen Lösungsmitteln. Halten Sie bei der Entsorgung der Abfälle ein geregeltes und genehmigtes Verfahren ein. Entsorgen Sie entzündliche, toxische und/oder infektiöse Substanzen keinesfalls über die öffentliche Kanalisation.

#### • Gefährliche Gase

Stellen Sie das HPLC-System in einem gut belüfteten Labor auf. Wenn die mobile Phase oder Probe flüchtige oder entzündliche Lösungsmittel enthält, müssen Sie sicherstellen, dass diese nicht in Ihren Arbeitsbereich gelangen. Vermeiden Sie offenes Feuer und Funken, wenn die mobile Phase oder Probe flüchtige oder entzündliche Stoffe enthält.

#### • Elektrostatische Entladung

Elektrostatische Entladung kann zu Funkenbildung führen und eine Brandgefahr darstellen. Beachten Sie, dass sich fließende Lösungsmittel in Kapillaren selbsttätig aufladen können. Besonders stark kann dieser Effekt in isolierenden Kapillaren und bei nicht leitenden Lösungsmitteln (beispielsweise reines Acetonitril) auftreten.

Ergreifen Sie geeignete Maßnahmen, um elektrostatische Aufladungen im Bereich des HPLC-Systems zu verhindern. Sorgen Sie beispielsweise für eine ausreichende Luftfeuchtigkeit und Belüftung im Labor, tragen Sie antistatische Schutzkleidung, vermeiden Sie die Ansammlung von Luftblasen in Abfallleitungen und verwenden Sie geerdete Abfallbehälter. Verwenden Sie nur nicht-leitende Kapillaren, um Lösungsmittel in den Abfallbehälter zu leiten. Elektrisch leitende Kapillaren sollten grundsätzlich geerdet sein.

#### Selbstentzündung von Lösungsmitteln

Verwenden Sie keine Lösungsmittel, deren Selbstentzündungstemperatur unter 150 °C liegt. Bei einer Undichtigkeit könnten sich diese Lösungsmittel an einer heißen Oberfläche selbst entzünden.

#### • Kapillaren, Kapillarverbindungen, offene Verbindungen

- Kapillaren, insbesondere nichtmetallische Kapillaren, können bersten, aus den Verschraubungen rutschen oder nicht eingeschraubt sein. Dies kann auch dazu führen, dass Substanzen aus den offenen Verbindungen spritzen.
- ◆ In einem UltiMate 3000-System kommen auch Komponenten aus PEEK<sup>™</sup> zum Einsatz. Dieses Polymer weist eine ausgezeichnete Beständigkeit gegen die meisten organischen Lösungsmittel auf. Es neigt jedoch dazu aufzuquellen, wenn es mit Trichlormethan (CHCl<sub>3</sub>), Dimethylsulfoxid (DMSO) oder Tetrahydrofuran (THF) in Kontakt kommt. Konzentrierte Säuren wie Schwefel- und Salpetersäure oder ein Gemisch aus Hexan, Ethylacetat und Methanol können PEEK angreifen. Beides kann dazu führen, dass Kapillaren undicht werden oder bersten. Die konzentrierten Säuren stellen bei kurzen Spülzyklen jedoch kein Problem dar.
- Verwenden Sie keine übermäßig beanspruchten, verbogenen, geknickten oder beschädigten Kapillaren.
- Kapillarverschraubungen können mit gefährlichen Substanzen kontaminiert sein oder es können gefährliche Substanzen an den offenen Verbindungen austreten.

- In einem UltiMate 3000 Bio RS-System kommen Viper-Systemkapillaren zum Einsatz, die aus der Nickel-Kobalt-Legierung MP35N<sup>®</sup> gefertigt sind. Hautkontakt mit diesem Material kann bei Personen, die gegen Nickel/Kobalt empfindlich sind, gegebenenfalls eine allergische Reaktion hervorrufen.
- Tragen Sie beim Umgang mit Fused Silica-Kapillaren immer eine Schutzbrille, z.B. bei der Installation oder zum Ablängen der Kapillaren.
- Ziehen Sie den Netzstecker, ehe Sie Abdeckungen am Gerät entfernen. Einige Bauteile im Innern des Gerätes können Spannung führen. Das Gehäuse darf nur vom Thermo Fisher Scientific-Kundendienst geöffnet werden.
- Tauschen Sie beschädigte Kommunikationskabel aus.
- Tauschen Sie beschädigte Netzkabel aus. Verwenden Sie nur die für das Gerät bereitgestellten Netzkabel.
- Ersetzen Sie durchgebrannte Sicherungen immer durch die von Thermo Fisher Scientific autorisierten Original-Ersatzsicherungen.
- Verwenden Sie ausschließlich die von Thermo Fisher Scientific für das Gerät autorisierten Original-Ersatz- und Zubehörteile.
- Schauen Sie niemals direkt in die LED für die Innenbeleuchtung des Autosamplers. Verwenden Sie keine lichtbündelnden Instrumente zur Betrachtung des Lichtstrahls. Die hohe Leuchtkraft der Lampe kann schädlich für die Augen sein
- Der Autosampler ist bei Auslieferung mit einer 2-Propanol-Lösung gefüllt. Achten Sie bei der ersten Inbetriebnahme darauf, dass das verwendete Lösungsmittel mit 2-Propanol mischbar ist. Ist dies nicht der Fall, stellen Sie das Lösungsmittel schrittweise um.
- Heben oder bewegen Sie den Autosampler nicht allein. Der Autosampler ist zu schwer und sperrig für eine Person allein.
- Wenn Sie den Autosampler anheben oder bewegen möchten, greifen Sie unter den Boden oder heben Sie das Gerät an den Seiten an. Schließen Sie ggf. vorher den Frontdeckel. Damit werden Schäden am Gerät vermieden.
- Der geöffnete Frontdeckel kann kein Gewicht aufnehmen. Legen Sie daher keine Gegenstände auf dem geöffneten Frontdeckel ab.
- Um Verletzungen zu vermeiden, greifen Sie während einer laufenden Analyse niemals in den Probenraum.
- Verwenden Sie nur handelsübliche Lösungsmittel mindestens in HPLC-Qualität und Puffer, die kompatibel mit den medienberührten Teilen des Autosamplers sind.
- Wenn eine Undichtigkeit auftritt, schalten Sie den Autosampler sofort aus und beheben Sie die Ursache für die Undichtigkeit.

- Beachten Sie bei längeren Betriebsunterbrechungen (= mehrere Tage) und wenn Sie den Autosampler für den Versand vorbereiten die Hinweise zur Außerbetriebnahme des Autosamplers (→ Seite 98).
- Setzen Sie den Autosampler nur entsprechend der bestimmungsgemäßen Nutzung des Gerätes und den Beschreibungen in dieser *Bedienungsanleitung* ein.
- Bewahren Sie die Bedienungsanleitung in Gerätenähe auf, so dass sie bei Bedarf schnell zur Hand ist.

### **1.3 Verwendungszweck des Autosamplers**

Das Gerät wurde ausschließlich für Forschungsaufgaben entwickelt. Es ist nicht für den Einsatz in diagnostischen Verfahren gedacht.

Es darf nur von qualifiziertem und berechtigtem Laborpersonal betrieben werden. Alle Benutzer müssen die Gefahren kennen, die vom Gerät und den verwendeten Substanzen ausgehen.

Der Autosampler wurde entwickelt für Laborforschungsaufgaben in der

- High-Performance Liquid Chromatography WPS-3000SL Analytical, WPS-3000SL Semiprep und WPS-3000TBSL Analytical
- Ultra-High Performance Liquid Chromatography WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS

Er ist Teil des UltiMate 3000-Systems, kann aber auch mit anderen Systemen betrieben werden, die über die entsprechenden Ein- und Ausgänge zur Steuerung des Moduls verfügen. Dazu wird ein Rechner mit USB-Port benötigt.

Der Autosampler wird über das Chromatographie-Management-System Chromeleon gesteuert. Als Teil des UltiMate 3000-Systems kann er auch mit anderen Datensystemen betrieben werden, wie

- Xcalibur<sup>™</sup>, Compass<sup>™</sup>/HyStar<sup>™</sup> oder Analyst<sup>®</sup>
   Dazu muss zusätzlich zum jeweiligen Datensystem die Software DCMS<sup>Link</sup> (Dionex Chromatography Mass Spectrometry Link) installiert werden.
- Empower<sup>TM</sup>

Dazu muss zusätzlich zum Datensystem die entsprechende Dionex Instrument Integration Software installiert werden.

Beachten Sie für die Verwendung des Autosampler Folgendes:

- Der Autosampler darf nur mit den von Thermo Fisher Scientific empfohlenen Zubehörund Ersatzteilen (→ Seite 173) und innerhalb seiner technischen Spezifikationen (→ Seite 163) betrieben werden.
- Verwenden Sie nur handelsübliche Lösungsmittel und Puffer, die kompatibel sind mit den medienberührten Teilen Ihres UltiMate 3000-Systems. Informationen zu den Materialien, aus denen die medienberührten Teile des Autosamplers gefertigt sind, finden Sie im Kapitel Technische Daten (→ Seite 163). Informationen zu den medienberührten Teilen der anderen Module Ihres UltiMate 3000-Systems finden Sie im Kapitel 'Technische Daten' in der Bedienungsanleitung für das jeweilige Modul.
- Achten Sie auf spezifische Eigenschaften der Lösungsmittel, wie Viskosität, Siedepunkt, UV-Absorption (UV/VIS-Detektor) und Brechungsindex (Brechungsindex-Detektor).

- pH-Bereich Abhängig vom Rotor Seal im Injektionsventil (→ Kap. 7.8.4, Seite 155).
- Pufferkonzentration
   WPS-3000SL Analytical, WPS-3000 Semiprep, WPS-3000TBSL Analytical,
   WPS-3000RS und WPS-3000TXRS:
   Typischerweise bis zu 1 mol/L (< 0,1 mol/L Chlorid-Ionen)
   WPS-3000TBRS:
   Typischerweise bis zu 1 mol/L ( ≤1 mol/L Chlorid-Ionen)
- Beachten Sie auch die Hinweise zur Lösungsmittelkompatibilität der anderen Module Ihres UltiMate 3000-Systems. Entsprechende Informationen finden Sie in den Bedienungsanleitungen zu den einzelnen Modulen.

Fragen zur bestimmungsgemäßen Nutzung des Gerätes beantworten wir Ihnen gern.

Für Schäden materieller oder immaterieller Art, die sich aus einer nicht bestimmungsgemäßen Verwendung bzw. unsachgemäßen Anwendung des Gerätes ergeben, kann Thermo Fisher Scientific keine Haftung übernehmen.

**Warnung:** Wenn das Gerät nicht entsprechend den Angaben von Thermo Fisher Scientific eingesetzt wird, kann der durch das Gerät bereitgestellte Schutz beeinträchtigt werden. Thermo Fisher Scientific übernimmt dann keine Verantwortung und haftet nicht für Verletzungen des Bedieners und/oder Schäden am Gerät. Wenn der Sicherheitsschutz des Gerätes zu irgendeinem Zeitpunkt nicht mehr gewährleistet ist, ist das Gerät von allen Stromquellen zu trennen und gegen jeden Betrieb zu sichern

STOP

# 2 Überblick

# 2.1 Kurzbeschreibung

Der Autosampler ist ein unentbehrlicher Baustein innerhalb des UltiMate 3000-Systems und bietet hohe Verfügbarkeit und Reproduzierbarkeit auch bei kleinsten Injektionsvolumina. Das Design wurde auf minimales Totvolumen und maximale Effizienz optimiert.

- Der Drehteller kann Probenfläschchen in verschiedenen Größen sowie Mikrotiter- und Deep-Well-Platten in einer beliebigen Kombination aufnehmen (→ Seite 23), die alle in derselben Autosampler-Sequenz verwendet werden können.
- Der Autosampler arbeitet nach dem "Split-Loop"-Injektionsprinzip, wodurch bei kurzer Injektionszeit (cycle time) ein variables Injektionsvolumen ermöglicht und ausgezeichnete Volumenreproduzierbarkeit sowie minimale Verschleppung gewährleistet werden (→ Seite 12).
- Zur Verringerung der Gradientenverzögerungszeit und Minimierung der Spülzeit unterstützt der Autosampler den Bypass-Modus (→ Seite 13).
- Zur Fehlererkennung stehen im Hinblick auf optimale Geräteleistung und Zuverlässigkeit diverse Sicherheits- und Überwachungsfunktionen zur Verfügung (→ Seite 29).
- Abhängig von den installierten Komponenten unterstützt der Autosampler Applikationen in der analytischen, Mikro- oder semipräparativen HPLC.
- Der Autosampler kann vollständig über das Chromatographie-Management-System Chromeleon gesteuert werden. Damit können Sie wahlfrei auf die Proben zugreifen, mehrere Standards in einer Probenserie verwenden sowie das Injektionsvolumen und die Anzahl der Replikate für jede Probe individuell festlegen.
- Jede Geräteversion steht auch mit integrierter Temperaturregelung zur Kühlung/Erwärmung der Proben zur Verfügung (→ Seite 16).
- Die fluidischen Komponenten im Autosampler sind leicht zugänglich, so dass Wartungsarbeiten schnell und zuverlässig durchgeführt werden können, während sich der Autosampler im UltiMate 3000-System befindet.
- Alle medienberührten Teile sind aus Materialien gefertigt, die eine ausgezeichnete Beständigkeit gegen die meisten eingesetzten Lösungsmittel und Puffer erreichen.
- Eine transparente Frontabdeckung (Best.-Nr. 6820.1427) ist als Option verfügbar. Damit können Sie den Probenraum des Autosamplers verschließen, zum Beispiel zum Schutz gegen Staub oder andere Partikel.

# 2.2 Funktionsprinzip

Die Abbildung illustriert das Funktionsprinzip des Autosamplers:



Abb. 1: Funktionsprinzip

Das Split-Loop-Injektionsprinzip beruht darauf, dass die Probennadel selbst Teil der Probenschleife ist: Während das Injektionsventil in der Position **Load** steht, wird die Probe mit Hilfe der Spritze durch die Probennadel in die Probenschleife gesaugt. Anschließend fährt die Probennadel in den Nadelsitz und das Injektionsventil wird in die Position **Inject** umgeschaltet. Das Laufmittel durchströmt nun die Probennadel und die Probenschleife und bringt die Probe auf die Trennsäule. Dadurch erübrigt sich das Spülen der Probennadel von innen. Außerdem wird eine exakte Dosierung variabler Probenvolumina ermöglicht. Dieses Prinzip ist verlustfrei. Die gesamte aufgezogene Probenmenge wird auch injiziert.

Der Autosampler enthält die folgenden fluidischen Komponenten:

Komponente	Beschreibung
Injektionsventil/ Probenschleife	Über das Injektionsventil und die Probenschleife wird die Probe in die mobile Phase eingebracht.
Probennadel	Die Probennadel nimmt die Probe auf und transportiert sie in die Probenschleife.
Waschport	Im Waschport kann das Äußere der Probennadel mit der Waschflüssigkeit gespült werden. Zusätzlich wird das Äußere der Probennadel nach dem Waschvorgang über die Membranpumpe mit Luft getrocknet.
	Wenn sich die Nadel in der untersten Position im Waschport befindet, wird die Nadelspitze direkt mit der Membranpumpe verbunden. Die Pumpe eines UltiMate 3000-Systems kann dann über die Membranpumpe entlüftet werden. Weitere Informationen dazu finden Sie im <i>Pumpenhandbuch</i> .

Komponente	Beschreibung
Nadelsitz	Der Nadelsitz dichtet die Nadelspitze gegen den Systemdruck ab. Eine spezielle Geometrie verhindert Verschleppung und dass Luft in den Hochdruckkreislauf injiziert wird.
Drehteller	Der Drehteller ist mit den Probenhaltern und/oder Mikrotiterplatten bestückt.
Spritze	Über die Spritze wird die Probe aus der Kavität bzw. dem Probenfläschchen in die Probenschleife gesaugt. Außerdem versorgt die Spritze den Waschport mit Reinigungslösung.
Pufferschleife	Die Pufferschleife schützt die Spritze vor Verunreinigungen durch die mobile Phase.
Waschflüssigkeit	Nach einer Injektion wird mit der Waschflüssigkeit die mobile Phase aus der Pufferschleife entfernt. Wurde eine Injektion unterbrochen, kann mit der Waschflüssigkeit die Probe aus der Probennadel und der Probenschleife entfernt werden.
Membranpumpe	Nach dem Waschvorgang wird das Äußere der Probennadel über die Membranpumpe mit Luft getrocknet. Wenn sich die Nadel in der untersten Position im Waschport befindet, wird
	die Nadelspitze direkt mit der Membranpumpe verbunden. Die Pumpe eines UltiMate 3000-Systems kann dann über die Membranpumpe entlüftet werden. Weitere Informationen dazu finden Sie im <i>Pumpenhandbuch</i> .
X und Z	Bezeichnen die möglichen Bewegungsrichtungen der Nadel.
Υ	Bezeichnet die mögliche Bewegungsrichtung des Drehtellers.

#### **Bypass-Modus**

Wie oben beschrieben, sind die Probennadel und die Probenschleife beim Split-Loop Injektionsprinzip Teil des chromatographischen Flusses, während sich das Injektionsventil in der Position **Inject** befindet (main pass). Das Laufmittel durchströmt die Probennadel und die Probenschleife und bringt die Probe auf die Trennsäule.

Um die Gradientenverzögerungszeit und die Spülzeit weiter verringern zu können, können Probennadel und Probenschleife aus dem chromatographischen Fluss herausgenommen werden. Dazu wird das Injektionsventil nach einer bestimmten Zeit in die Position **Load** geschaltet (= Bypass-Modus). Das Laufmittel bringt die Probe nun direkt auf die Trennsäule auf. Die Zeit, nach welcher das Injektionsventil umgeschaltet werden soll, können Sie in Chromeleon über den **FlushOutFactor** festlegen. Geben Sie dort das n-fache des Injektionsvolumens an. Das "Flush out"-Volumen ist das Volumen, das benötigt wird, um die Probe auf die Trennsäule aufzubringen, wenn sich das Injektionsventil in der Position **Inject** befindet. Es hängt vom Injektionsvolumen und dem Volumen aller Kapillaren ab, die die Probe durchströmen muss, ehe sie das Injektionsventil verlässt. Chromeleon rechnet diesen Faktor unter Berücksichtigung der Flussrate in die Zeit um, nach der das Injektionsventil in die Position **Load** geschaltet wird.

Außerdem kann im Bypass-Modus der Probendurchsatz erhöht werden, da bereits vor dem Ende der Analyse der aktuellen Probe mit der Vorbereitung der nächsten Probe begonnen werden kann. Weitere Informationen finden Sie in der *Chromeleon-Hilfe*.

# 2.3 Autosamplerkonfigurationen

#### 2.3.1 Übersicht

Der Autosampler ist in den in der Tabelle genannten Konfigurationen verfügbar. Informationen zu den optional erhältlichen Komponenten finden Sie auf Seite 178.

WPS-3000SL	BestNr.
<ul> <li>WPS-3000SL Analytical (ohne Probenthermostatisierung) mit</li> <li>100-μL-Spritze</li> <li>100-μL-Probenschleife</li> <li>100-μL-Pufferschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet f ür einen Druck &lt; 62 MPa (9000 psi)</li> </ul>	5822.0010
WPS-3000TSL Analytical, wie WPS-3000SL Analytical, jedoch mit Probenthermostatisierung	5822.0020
<ul> <li>WPS-3000SL Semiprep (ohne Probenthermostatisierung) mit</li> <li>2500-μL-Spritze</li> <li>2500-μL-Probenschleife</li> <li>&gt;250-μL-Pufferschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet f ür einen Druck &lt; 62 MPa (9000 psi)</li> </ul>	5822.0018
WPS-3000TSL Semiprep wie WPS-3000SL Semiprep, jedoch mit Probenthermostatisierung	5822.0028

WPS-3000TBSL	BestNr.
<ul> <li>WPS-3000TBSL Analytical (für elektrochemische Detektion) mit</li> <li>100-μL-Spritze</li> <li>100-μL-Probenschleife</li> <li>100-μL-Pufferschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet für einen Druck &lt; 34 MPa (5000 psi)</li> <li>Probenthermostatisierung</li> </ul>	

WPS-3000RS	BestNr.
<ul> <li>WPS-3000RS (ohne Probenthermostatisierung) mit</li> <li>100-μL-Spritze</li> <li>100-μL-Probenschleife</li> <li>100-μL-Pufferschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet f ür einen Druck &lt; 103 MPa (15000 psi)</li> </ul>	5840.0010
WPS-3000TRS, wie WPS-3000RS, jedoch mit Probenthermostatisierung	5840.0020
Der WPS-3000(T)RS kann für Anwendungen in der Mikro-HPLC optimiert werden	

 $(\rightarrow$  Seite 161).

WPS-3000TBRS	BestNr.
<ul> <li>WPS-3000TBRS (biokompatibler Autosampler) mit</li> <li>100-μL-Spritze</li> <li>25-μL-Probenschleife</li> <li>100-μL-Pufferschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet f ür einen Druck &lt; 103 MPa (15000 psi)</li> <li>Probenthermostatisierung</li> </ul>	5841.0020

WPS-3000TXRS	BestNr.
WPS-3000TXRS mit	5843.0020
- 25-µL-Probenschleife	
- Injektionsventil (2 Positionen, 6 Ports)	
geeignet für einen Druck < 125 MPa (18130 psi) - Probenthermostatisierung	

#### 2.3.2 Autosampler mit Probenthermostatisierung

Autosampler mit Probenthermostatisierung (angezeigt durch das 'T' im Namen) verfügen über thermoelektrische (Peltier-) Elemente, die den Drehteller und die darin befindlichen Komponenten kühlen oder erwärmen und so erlauben eine präzise Temperierung der Proben erlauben.

Ausgehend von der Umgebungstemperatur können die Proben um max. 22 °C gekühlt werden. Die untere Temperaturgrenze für die Kühlung beträgt +4 °C. Darüber hinaus können die Proben auf max. +45 °C erwärmt werden.

Zum Betrieb sind keine weiteren Anschlüsse wie Kühlwasser, Druckluft oder ähnliches erforderlich.

Beachten Sie die folgenden Hinweise:

- Zur Ableitung von eventuell entstehendem Kondenswasser können Sie rechts hinten auf der Unterseite des Gerätes einen Ablaufschlauch anschließen (→ Abb. 18, Seite 54). Der Autosampler verfügt über eine aktive Kondenswasserdrainage (über eine Pumpe).
- Achten Sie beim Anschluss des Ablaufschlauches darauf, dass dieser an keiner Stelle höher liegt als der Anschluss-Stutzen. Bei Missachtung kann es zu einer Beschädigung des Gerätes kommen, da Kondenswasser in das Gerät fließen kann.
- Wischen Sie bei starker Kondenswasserbildung den Drehteller von Zeit zu Zeit aus, um Verunreinigungen zum vermeiden.

#### 2.3.3 Kombinationsmöglichkeiten von Spritzen und Schleifen

Die nachfolgende Tabelle gibt einen kurzen Überblick über die empfohlenen Kombinationen von Spritze, Pufferschleife und Probenschleife.

<b>Spritze</b> (BestNr. siehe Seite 121)	<b>Pufferschleife</b> (BestNr. siehe Seite 124)	<b>Probenschleife<sup>1</sup></b> (BestNr. siehe Seite 131)
25 μL	100 µL	25 μL (V = 40 μL)
100 µL	100 μL	25 $\mu$ L (V = 40 $\mu$ L) oder 100 $\mu$ L (V = 130 $\mu$ L bzw. V = 150 $\mu$ L, je nach Probenschleifentyp)
250 μL	>250 µL	250 μL (V = 344 μL)
500 µL	>250 μL	500 μL (V = 667 μL) 1000 μL <sup>2</sup> (V = 1618 μL)
1000 µL	>250 μL	$1000 \ \mu L^2 \ (V = 1618 \ \mu L)$
2500 μL	>250 µL	$2500 \ \mu L^2 \ (V = 3300 \ \mu L)$

<sup>1</sup> Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen.

<sup>2</sup> Diese Probenschleifen sind für den WPS-3000SL Semiprep gedacht. Wenn Sie diese Probenschleifen mit einem WPS-3000SL Analytical verwenden möchten, ist ein Upgrade auf die semipräparative Autosamplerkonfiguration erforderlich. Wenden Sie sich in diesem Fall an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.

Einige in der Tabelle der genannten Kombinationen stehen auch als Injektionsvolumen-Kits zur Verfügung (→ Seite 178).

# 2.4 Innenansicht



Abb. 2: Innenansicht

Nr.	Beschreibung		
1	Frontdeckel (nach oben aufgeklappt)		
2	Spritzenventil ( $\rightarrow$ Seite 53)		
3	Beleuchtung des Probenraums, hier durch Frontdeckel verdeckt ( $\rightarrow$ Seite 25)		
4	Schlauchverbindung Spritzenventil - Waschport (Versorgung Waschport)		
5	Spritze		
6	Position für optionale Waschflasche (für 125 mL Waschflüssigkeit; → Seite 50)		
7	Menüstift		
8	Deckel Drehteller Sie können auf die Proben im Drehteller zugreifen, wenn Sie den Deckel am schwarzen Griff nach links schwenken. Um den Drehteller zu verschließen, schwenken Sie den Deckel ganz nach rechts, bis Kontakt zum Nadelarm besteht. Zum Aufziehen der Probe wird der Deckel durch die Bewegungen des Nadelarms automatisch geöffnet und geschlossen.		
9	Pufferschleife		
10	Leaksensor		
11	Auffangwanne ( $\rightarrow$ Detailansicht in Abb. 3, Seite 19)		
12	Injektionsventil ( $\rightarrow$ Seite 27)		
13	Probenschleife		



Abb. 3: Auffangwanne

Nr.	Beschreibung	
1	Waschport	
2	Injektionsventil ( $\rightarrow$ Seite 27)	
3	Pufferschleife (Port 1 am Injektionsventil)	
4	Probenschleife (Port 6 am Injektionsventil)	
5	Kapillare vom Injektionsventil (Port 2) zum Waschport (Injektor-Waste-Kapillare)	
6	Waschport (Unterteil)	
7	Kapillare zur Ableitung der Flüssigkeiten aus dem Waschport (Drainage)	
8	Schlauchverbindung Spritzenventil - Waschport (Versorgung Waschport)	
9	Durchlass zum Säulenthermostaten Führen Sie die Kapillare von Injektionsventil zum Eingang der Säule durch diesen Durchlass. Dies ist der kürzeste Weg zum Säulenthermostaten.	
10	Leaksensor ( $\rightarrow$ Seite 26)	
11	Kapillare vom Nadelsitz zum Injektionsventil (Port 3) (Nadelsitzkapillare)	
12	Nadelsitz	
13	Niederhalter mit Sensor zur automatischen Probenfläschchen- bzw. Mikrotiterplatten- und Deep-Well-Plate-Erkennung (→ Seite 95).	
14	Probennadel ( $\rightarrow$ Seite 125)	

### 2.5 Gerätevorderseite



Abb. 4: Gerätevorderseite

Nr.	Element	Funktion		
1	Display	<ul> <li>Zeigt verschiedene Informationen zum Autosampler an, z.B.:</li> <li>allgemeine Informationen beim Einschalten des Autosamplers (→ Seite 63)</li> <li>Statusanzeige (→ Seite 64)</li> <li>verschiedene Funktionen und Menüs (→ Seiten 73 und 75)</li> <li>Meldungen (→ Seite 104)</li> </ul>		
2	Standby-Taste	schaltet den Autosampler in den Standby-Modus (die LED leuchtet). Erneutes Drücken der Standby-Taste hebt den Standby-Modus auf (die LED leuchtet nicht). <b>Hinweis</b> : Damit der Autosampler den Modus ändert, muss die Standby- Taste ca.1 Sekunde lang gedrückt werden.		
3	Status-LEDs	-LEDs		
	Power	Die LED leuchtet blau, wenn der Autosampler eingeschaltet ist.		
	Connected	Die LED leuchtet grün, wenn der Autosampler mit Chromeleon verbunden ist.		
	Status	Die LED leuchtet standardmäßig grün; der Autosampler ist betriebsbereit. Die LED leuchtet rot, wenn ein Fehler erkannt wurde. Autosampler mit Probenthermostatisierung: Die LED leuchtet gelb, wenn die Solltemperatur noch nicht erreicht ist.		

# 2.6 Geräterückseite



Abb. 5: Anschlüsse auf der Geräterückseite (hier Autosampler mit Probenthermostatisierung)

Nr.	Beschreibung
1	Netzschalter ( $\rightarrow$ Seite 22)
2	Sicherungsschlitten ( $\rightarrow$ Seite 22)
3	Netzanschluss ( $\rightarrow$ Seite 36)
4	Typenschild
5	USB 1.1-Schnittstelle (Universal Serial Bus): Anschluss des Autosamplers an den Chromeleon-Rechner ( $\rightarrow$ Seite 22)
6	Interner USB-Hub (3 USB 1.1 Anschlussports) Abhängig von der Konfiguration des UltiMate 3000-Systems, gegebenenfalls Anschluss je eines weiteren Moduls der UltiMate 3000-Serie oder Anschluss eines USB-Hubs (→ Seite 22)
7	Digital I/O zur Kommunikation mit externen Geräten unter Chromeleon ( $\rightarrow$ Seite 22)
8	(Nur Autosampler mit Probenthermostatisierung) Ventilatoren

#### 2.6.1 Netzschalter

Der Hauptnetzschalter befindet sich auf der Geräterückseite. Über diesen Schalter wird der Autosamplerein- und ausgeschaltet.

#### 2.6.2 Sicherungsschlitten

Im Sicherungsschlitten befinden sich zwei Sicherungen. Abhängig vom Autosamplertyp sind dies zwei Sicherungen mit 2 A, 250 V (Autosampler ohne Probenthermostatisierung) oder 4 A, 250 V (Autosampler mit Probenthermostatisierung). Informationen zum Sicherungswechsel finden Sie auf Seite 158.

#### 2.6.3 USB-Anschluss (USB 1.1)

Der Autosampler kann über eine USB-Verbindung (USB = Universal Serial Bus) über das Chromatographie-Management-System Chromeleon gesteuert werden. Die Daten werden digital über das entsprechende USB-Anschlusskabel übertragen ( $\rightarrow$  Seite 35).

Je nach Konfiguration des UltiMate 3000-Systems, können Sie über den internen 3-Port-USB-Hub ( $\rightarrow$  Abb. 5, Nr. 6) weitere Module der UltiMate 3000-Reihe oder externe USB-Hubs anschließen. Der Anschluss von Geräten mit USB 2.0 ist nicht möglich.

**Vorsicht:** Thermo Fisher Scientific empfiehlt, die USB-Anschlüsse nur für den Anschluss von Dionex-Geräten zu verwenden. Werden Geräte anderer Anbieter angeschlossen, kann Thermo Fisher Scientific keine Funktionsgarantie übernehmen.

Informationen zum Anschluss des Autosamplers an den Chromeleon-Rechner finden Sie in den Kapiteln 3.4.1 und 3.4.2 ( $\rightarrow$  Seite 35).

#### 2.6.4 Digital I/O

Die 6-poligen Mini-DIN-Buchsen können unter Chromeleon zur Übertragung von digitalen Signalen mit externen Geräten verwendet werden. Die Übertragung der Signale erfolgt über ein entsprechendes Signalkabel (Best.-Nr. 6000.1004). Informationen zum Kabel und zum Anschluss finden Sie auf Seite 36. Informationen zur Anschlussbelegung der Buchsen und des Kabels finden Sie auf Seite 191.

Wenn das UltiMate 3000-System eine Pumpe LPG-3400XRS enthält, muss die Pumpe mit einem Digital I/O-Port des Autosamplers verbunden werden, wenn der Injektionsbefehl des Autosamplers mit den Kolbenhüben der Pumpe synchronisiert werden soll. Einzelheiten finden Sie in der *Bedienungsanleitung* zur Pumpe.

### 2.7 Drehteller, Probenhalter und Mikrotiterplatten

Der Drehteller ist in drei Segmente (rot, grün und blau) unterteilt und kann drei Probenhalter (Trays) in beliebiger Kombination aufnehmen. Informationen zur Installation der Probenhalter finden Sie auf Seite 57.



Abb. 6: Probenhalter (Beispiele)

Folgende Probenhalter stehen für Probenfläschchen zur Verfügung:

Beschreibung	BestNr.
Probenhalter für 40 Fläschchen (1,8 mL oder 2 mL)	6820.4070
Probenhalter für 40 Fläschchen (konisch; 1,1 mL)	6820.4087
Probenhalter für 40 Fläschchen (Eppendorf, 1,5 mL)	6820.4094
Probenhalter für 40 Fläschchen (Eppendorf, 0,5 mL)	6820.4096
Probenhalter für 10 Fläschchen (10 mL)	6820.4086
Probenhalter für 22 Fläschchen (4 mL)	6820.4084
Probenhalter für 72 Fläschchen (zylindrisch, 1,2 mL)	6820.4090
Probenhalter für72 Fläschchen (zylindrisch, 0,3 mL)	6820.4091
Probenhalter für 72 Mikrodialyse-Fläschchen (0,3 mL)	6820.4097

Beschreibung	BestNr.
Probenhalter für Standard-Mikrotiterplatte (Platten-Höhe: 12 - 24 mm)	6820.4070, 6820.4084, 6820.4086, 6820.4087, 6820.4090 oder 6820.4091
Probenhalter für Deep-Well-Plate (Plate-Höhe: 34 - 46 mm)	6820.4079
Probenhalter für Deep-Well-Plate (Plate-Höhe: 30 - 36 mm)	6820.4083
Probenhalter für Deep-Well-Plate (Plate-Höhe: 20 – 32 mm)	6820.4089
Probenhalter für Low-Well-PCR-Plates (Plate-Höhe: 8 - 12 mm)	6820.4070, 6820.4086, 6820.4087, 6820.4090 <i>oder</i> 6820.4091
	Zusätzlich ist ein Adapter erforderlich (BestNr. 6820.4088).

Folgende Probenhalter stehen für Mikrotiterplatten zur Verfügung:

Darüber hinaus sind an drei Positionen im Drehteller Halter für jeweils 5 Probenfläschchen à 10 mL, z.B. zur Aufnahme der Reagensflüssigkeiten, installiert ( $\rightarrow$  Seite 59).

Probengefäß (Min Max. Höhe)	Beschreibung	BestNr.
Probenfläschchen (31 - 43 mm)	Zylindrisches Probenfläschchen bis 1,8 mL (A.D. < 12 mm), z.B. Probenfläschchen 250 µL Kappe und Septum für 250-µL-Probenfläschchen für 40-Positionen-Probenhalter (BestNr. 6820.4070)	6820.0029 6820.0028
1,2-mL-Probenfläschchen (31 - 43 mm)	Zylindrisches Probenfläschchen Crimpkappe Septum (Silikon/PTFE, geschlitzt) für 72-Positionen-Probenhalter (BestNr. 6820.4090)	6000.0062 6000.0064 6000.0061
1,1-mL-Probenfläschchen (31-43 mm)	Konisches Probenfläschchen Crimpkappe und geschlitztem Silikon/PTFE Septum für 40-Positionen-Probenhalter (BestNr. 6820.4087)	6000.0077 6000.0076
1,8-mL-Probenfläschchen (31 - 43 mm)	Probenfläschchen Crimpkappe und Septum für 40-Positionen-Probenhalter (BestNr. 6820.4070)	6000.0072 6000.0071
2-mL-Probenfläschchen (31 - 43 mm)	Zylindrisches Probenfläschchen Schraubkappe und Silikondichtung für 40-Positionen-Probenhalter (BestNr. 6820.4070)	6000.0060 6000.0057 6000.0058

Informationen zu den von Thermo Fisher Scientific empfohlenen Mikrotiterplatten und Probenfläschchen sowie zu den erforderlichen Gefäßhöhen finden Sie in der Tabelle:

Probengefäß (Min Max. Höhe)	Beschreibung	BestNr.
4-mL-Probenfläschchen (36 - 48 mm)	Zylindrisches Probenfläschchen Schraubkappe und Silikon/PTFE-Septum für 22-Positionen Probenhalter (BestNr. 6820.4084)	6000.0074 6000.0073 6000.0075
10-mL-Probenfläschchen (38 - 50 mm)	Zylindrisches Probenfläschchen mit Kappe und Septum für 5-Positionen Halter (BestNr. 6820.4073)	6820.0023
Standard-Mikrotiterplatte (12 - 24 mm)	96 Kavitäten 384 Kavitäten Setzen Sie diese Platten auf einen der auf Seite 23 genannten Probenhalter auf.	6820.4100 6820.4110
Deep-Well-Plate (34 - 46 mm)	<ul> <li>96 Kavitäten</li> <li>384 Kavitäten</li> <li>Für die Deep-Well-Plates wird ein spezieller</li> <li>Plattenhalter benötigt (BestNr. 6820.4079).</li> <li>Installieren Sie Deep-Well-Plates <i>nicht</i> auf einem der anderen Probenhalter.</li> <li>Es können auch Deep-Well-Plates mit 24 Kavitäten verwendet werden. Dazu müssen Sie den Plattenhalter (BestNr. 6820.4079) <i>und</i> den Niederhalter-Adapter (BestNr. 6820.2402) installieren <i>und</i> den automatischen Tray-Test (→ Seite 94) ausschalten.</li> </ul>	6820.4101 Auf Anfrage

Beachten Sie auch Folgendes:

- Eine Verwendung anderer Septen kann zum Verschluss von Kapillaren im Gerät führen oder die chromatographischen Ergebnisse beeinflussen.
- Eppendorf-Fläschchen müssen offen sein.
- Bei Verwendung von 10-mL-Probenfläschchen *ohne* Kappe wird die Installation des Niederhalter-Adapters (Best.-Nr. 6820.2402) empfohlen. Wenn Sie den Adapter installieren, *müssen* Sie den automatischen Tray-Test ausschalten (→ Seite 94).
- Wenn Sie andere als die oben genannten Probengefäße verwenden möchten, beachten Sie die in der Tabelle genannten erforderlichen Gefäßhöhen.

### 2.8 Probenraumbeleuchtung

An der Innenseite des Frontdeckels sind zwei Lampen installiert, die den Probenraum während des Autosampler-Betriebs beleuchten. Um bei einem Eingriff in den Probenraum Verletzungen zu vermeiden, signalisieren sie darüber hinaus eine Bewegung des Nadelarms. Die Lampen schalten kurzzeitig aus und wieder ein, ehe der Nadelarm in eine andere Position fährt.

Die Probenraumbeleuchtung ist standardmäßig aktiviert. Wenn Ihre Proben lichtempfindlich sind, setzen Sie das Property **TrayIllumination** in Chromeleon auf **Off**.

### 2.9 Fluidische Anschlüsse

Die fluidischen Anschlüsse im Autosampler sind leicht zugänglich ( $\rightarrow$  Abb. 2, Seite 18). Klappen Sie den Frontdeckel nach oben auf. Der Deckel arretiert in der obersten Position.

**Vorsicht:** Der geöffnete Frontdeckel kann kein Gewicht aufnehmen. Legen Sie daher keine Gegenstände auf dem geöffneten Frontdeckel ab.

Verwenden Sie für die fluidischen Anschlüsse nur die mitgelieferten, fertig konfektionierten Anschlusskapillaren bzw. Original-Dionex-Ersatzkapillaren. Verwenden Sie benutzte Fittingverbindungen nur für dieselbe Kapillarverbindung wieder, um ein erhöhtes Totvolumen zu vermeiden.

### 2.10 Leaksensor

Der Autosampler verfügt über einen Leaksensor, über den Undichtigkeiten im Autosampler erkannt werden. Wenn die Leakerkennung eingeschaltet ist, spricht der Leaksensor an, wenn sich Flüssigkeit in der Auffangwanne unter den fluidischen Verbindungen sammelt. Die LED **Status** auf der Gerätevorderseite leuchtet rot. Es erscheint eine Meldung auf dem Gerätedisplay und im Chromeleon Audit Trail. Je nach Einstellung ertönt außerdem ein akustisches Signal ( $\rightarrow$  Seite 86).

Wenn der Leaksensor angesprochen hat, beseitigen Sie die Undichtigkeit und trocknen Sie den Leaksensor ( $\rightarrow$  Seite 120). Die LED **Status** bleibt rot, bis der Sensor trocken ist. Um die Meldung auf dem Gerätedisplay zu löschen, wählen Sie **Clear** auf der Navigationsleiste ( $\rightarrow$  Seite 77).

# 2.11 Injektionsventil

Der Autosampler ist mit einem Injektionsventil (2 Positionen, 6 Ports) ausgestattet, das für Anwendungen mit folgendem Druck geeignet ist:

Autosampler	Druck
WPS-3000SL (Analytical und Semiprep)	< 62 MPa (9000 psi)
WPS-3000TBSL Analytical	< 34 MPa (5000 psi)
WPS-3000RS	< 103 MPa (15000 psi)
WPS-3000TBRS	< 103 MPa (15000 psi)
WPS-3000TXRS	< 125 MPa (18130 psi)

Das Ventil hat zwei Betriebspositionen: Load und Inject. In der Position Load wird die Probe über die Probennadel in die Probenschleife gesaugt. Der Anschluss für die Pumpe ist direkt mit dem Anschluss für die Säule verbunden (Bypass-Modus). Wenn das Injektionsventil in die Position Inject geschaltet wird, wird die Probe in den Hochdruckkreislauf zur Analyse auf die Trennsäule aufgebracht.

Die Komponenten sind wie folgt am Ventil angeschlossen:

Port	Anschluss von	Port	Anschluss von
1	Spritze/Pufferschleife	4	Kapillare zur Säule
2	Injektor-Waste-Kapillare	5	Kapillare von der Pumpe
3	Nadelsitzkapillare	6	Probennadel/Probenschleife

Informationen zur Wartung des Ventils finden Sie im Kapitel 7.8 ( $\rightarrow$  Seite 148).

# 2.12 Steuerung über Chromeleon

Der Autosampler kann über das Chromatographie-Management-System Chromeleon gesteuert werden. Voraussetzung hierfür ist eine geeignete Chromeleon-Version sowie eine Lizenz der Klasse **Timebase Class 1**.

Der Autosampler kann auf zweierlei Art über Chromeleon gesteuert werden:

• Direkt

Beim direkten Betrieb wählen Sie die Parameter und Befehle im Dialogfenster **Commands** (F8-Box) aus. Direkte Befehle werden mit der Eingabe ausgeführt. Für den Routinebetrieb stehen die meisten Parameter und Befehle auch in einem Steuerfenster zur Verfügung. Weitere Informationen zum direkten Betrieb finden Sie auf Seite 68.

• Automatisch

Beim automatischen Betrieb erstellen Sie ein Steuerprogramm (PGM-File). Dabei handelt es sich um eine Liste von Steuerbefehlen, die in zeitlicher Abfolge ausgeführt werden. Über diese Liste wird der Autosampler automatisch gesteuert. Ein Programm können Sie automatisch mit einem Software-Assistenten erstellen oder manuell, indem Sie ein vorhandenes Programm editieren. Weitere Informationen zum automatischen Betrieb finden Sie auf Seite 71.

**Hinweis:** Alle Software-Beschreibungen in dieser Bedienungsanleitung beziehen sich auf *Chromeleon 6.80*.

Wenn Sie den Autosampler mit *Chromeleon 7* betreiben möchten, finden Sie Informationen zu den entsprechenden Arbeitsabläufen in den folgenden Dokumenten (alle im Lieferumfang von Chromeleon 7 enthalten):

- *Online-Hilfe zu Chromeleon* 7—bietet umfangreiche Informationen und ausführliches Referenzmaterial zu allen Aspekten der Software.
- *Quick Start Guide*—beschreibt die wichtigsten Elemente der Benutzeroberfläche und führt Sie schrittweise durch die wichtigsten Arbeitsabläufe.
- *Referenzkarte*—beschreibt die wichtigsten Arbeitsabläufe in Kurzform.
- *Installation Guide*—bietet grundlegende Informationen zur Installation und Konfiguration von Geräten. Spezifische Informationen zu einzelnen Geräten finden Sie in der Hilfe zum *Chromeleon 7 Instrument Configuration Manager.*

Beachten Sie auch, dass

- Chromeleon 7 eine andere Terminologie verwendet als Chromeleon 6.80. Informationen hierzu finden Sie in dem Dokument 'Glossary -Chromeleon 7' (enthalten im Ordner 'Documents' der Chromeleon 7-Installation).
- Gegebenenfalls einige Funktionen in Chromeleon 7 noch nicht unterstützt werden.
## 2.13 Wellness, Predictive Performance und Diagnose

Wellness überwacht das "Wohlbefinden" des Autosamplers. Damit aus kleinen Problemen keine großen Probleme werden, verfügt der Autosampler über verschiedene Sicherheitsfunktionen zur Überprüfung seiner Leistung und Zuverlässigkeit. Dazu gehören:

- Interne Überwachung aller mechanischen Operationen
- Automatischer Selbsttest beim Einschalten
- Automatische Mikrotiterplatten- und Probenfläschchenerkennung ( $\rightarrow$  Seite 95)
- Leaksensor ( $\rightarrow$  Seite 26)
- Überwachung der Kühlfunktion durch drei Temperaturfühler

Wird ein Fehler gefunden, leuchtet die LED **Status** auf der Gerätevorderseite rot und eine Meldung blinkt auf dem Gerätedisplay ( $\rightarrow$  Seite 104).

Wird der Autosampler über Chromeleon betrieben, stehen zusätzlich Funktionen zur aktiven Überwachung von Verschleißteilen (= Predictive Performance;  $\rightarrow$  Seite 88) sowie verschiedene Diagnosetests ( $\rightarrow$  Seite 89) zur Verfügung, mit denen der Autosampler und verschiedene seiner Komponenten auf ihre Funktionsfähigkeit hin überprüft werden können.

# 3 Installation

## 3.1 Anforderungen an den Standort

Der Standort muss die folgenden Anforderungen erfüllen:

- Der Hauptnetzschalter und der Netzanschluss befinden sich auf der Geräterückseite. Stellen Sie sicher, dass
  - der Hauptnetzschalter jederzeit einfach und frei zugänglich ist.
  - das Netzkabel des Gerätes einfach zugänglich ist und jederzeit vom Stromnetz getrennt werden kann. Lassen Sie hinter dem Gerät ausreichend Platz, damit das Netzkabel herausgezogen werden kann.
- Der Standort muss die in den technischen Daten (→ Seite 163) unter Leistungsaufnahme und Umgebungsbedingungen genannten Spezifikationen erfüllen.
- Stellen Sie den Autosampler auf eine stabile und vibrationsfreie Unterlage.
- Der Untergrund muss lösungsmittelresistent sein.
- Die Umgebungstemperatur sollte möglichst konstant sein.
- Vermeiden Sie direkte Sonneneinstrahlung und hohe Luftfeuchtigkeit.
- Achten Sie bei der Aufstellung des Autosamplers darauf, dass der Platz hinter und neben dem Gerät ausreichend für die Belüftung ist.

## 3.2 Auspacken

Alle Autosampler werden vor dem Versand sowohl elektrisch als auch mechanisch sorgfältig geprüft. Nach dem Auspacken prüfen Sie den Lieferumfang auf offensichtliche Anzeichen mechanischer Beschädigungen, die auf dem Transportweg aufgetreten sein könnten.

- **Hinweise:** 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. Bewahren Sie die Originalverpackung auf. Sie ist die bestmögliche Verpackung für den Transport des Gerätes (z.B. im Reparaturfall). Eine Gerätegarantie wird nur übernommen, wenn das Gerät in der Originalverpackung eingeschickt wird.
- 1. Stellen Sie den Karton auf den Boden und entnehmen Sie den Zubehörkarton sowie das Netzkabel. Die beweglichen Teile für den Drehteller (Probenhalter, Mikrotiterplatten) sind separat verpackt.
- 2. Heben Sie das Gerät vorsichtig aus dem Karton, und stellen Sie es auf eine stabile Unterlage.
  - ▲ Vorsicht: Heben oder bewegen Sie den Autosampler nicht allein. Der Autosampler ist zu schwer und sperrig für eine Person allein. Um ein Herunterfallen des Gerätes zu vermeiden, greifen Sie den Autosampler an der Geräteunterseite und heben Sie ihn zusammen mit den seitlichen Verpackungsteilen aus dem Karton. Wenn Sie den Autosampler anheben oder bewegen möchten, greifen Sie immer unter den Boden oder heben Sie das Gerät an den Seiten an.
- 3. Entfernen Sie die Verpackungsteile und die Kunststoff-Folie.
- 4. Entfernen Sie die Verpackungsschaumteile, die den Autosampler während des Transports schützen:
  - Entnehmen Sie das Schaumteil, das den Nadelarm und Drehteller sichert.
  - Zwei Schaumteile arretieren den Frontdeckel. Klappen Sie den Frontdeckel nach oben auf und entnehmen Sie die beiden Schaumteile.
- 5. Während des Transports kann sich durch große Temperaturunterschiede Kondenswasser im Gerät bilden. Temperieren Sie daher den Autosampler nach dem Auspacken mindestens 4 Stunden, damit das Kondenswasser entfernt wird. Schließen Sie den Autosampler dabei nicht an die Stromversorgung an. Bestehen offensichtliche Zweifel an der vollständigen Auflösung des Kondenswassers, ist der Autosampler so lange im elektrisch nicht angeschlossenen Zustand zu akklimatisieren, bis sich Niederschläge vollständig verflüchtigt haben.

## 3.3 Position des Autosamplers im UltiMate 3000-System

Wenn Sie den Autosampler innerhalb eines UltiMate 3000-Systems betreiben, stellen Sie die Module des Systems übereinander ( $\rightarrow$  Abb. 7) und verbinden Sie diese auf der Rückseite miteinander ( $\rightarrow$  Abb. 8). Der individuelle Systemaufbau hängt von der Anwendung ab.



Abb. 7: Beispiel für die Aufstellung eines UltiMate 3000-Systems



Abb. 8: Beispiel für die Verbindung der Module auf der Rückseite eines UltiMate 3000-Systems

Alle Module eines UltiMate 3000-Systems (mit Ausnahme des Solvent Racks) können auch direkt über die USB-Schnittstelle mit dem Chromeleon-Rechner verbunden werden. Thermo Fisher Scientific empfiehlt jedoch, alle Module untereinander zu verbinden und nur eine Verbindung zum Rechner zu führen.

Bei Systemen mit DAD-3000(RS) oder MWD-3000(RS) ist dies *nur* vom Detektor aus möglich.

Bei Systemen mit VWD 3x00(RS) ist dies nur von der Pumpe aus möglich.

## 3.4 Verbinden des Autosamplers

#### 3.4.1 Allgemeine Informationen

Wenn Sie den Autosampler mit Chromeleon betreiben möchten

*Bevor* Sie den Autosampler mit dem Chromeleon-Rechner verbinden und den Autosampler einschalten, vergewissern Sie sich, dass die Chromeleon-Software auf dem Rechner installiert und der Lizenzcode eingegeben ist. Nur wenn die Chromeleon-Software zuerst installiert ist und der Autosampler danach mit dem Rechner verbunden und eingeschaltet wird, wird der USB-Treiber für den Autosampler automatisch geladen. Nur dann kann das Windows<sup>®</sup>-Betriebssystem den Autosampler erkennen, wenn dieser eingeschaltet wird.

#### 3.4.2 Anschluss des USB-Kabels

Verbinden Sie den Autosampler über die USB-Anschlüsse ( $\rightarrow$  Abb. 5, Seite 21) auf der Geräterückseite mit dem Chromeleon-Rechner. Wählen Sie eine der folgenden Alternativen:

- Verbinden Sie den Autosampler direkt mit dem USB-Anschluss des Chromeleon-Rechners.
- Verbinden Sie den Autosampler mit dem internen USB-Anschluss eines anderen Moduls innerhalb des UltiMate 3000-Systems, das mit dem Chromeleon-Rechner verbunden ist (→ Seite 34). Der Anschluss von Geräten mit USB 2.0 ist nicht möglich.
- Verbinden Sie den Autosampler über einen externen USB-Hub mit dem Chromeleon-Rechner.

Um eine einwandfreie Funktion der Verbindung zu gewährleisten, verwenden Sie nur diese Kabel für die Verbindung:

USB-Kabel	BestNr.
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 1 m)	6035.9035 (im Autosampler-Zubehör enthalten)
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 5 m)	6911.0002

# **i** Hinweis:

Die Länge der USB-Verbindung zum Rechner oder nächsten USB-Hub darf 5 m nicht überschreiten.

#### 3.4.3 Anschluss des Netzkabels

Verbinden Sie die Netzbuchse auf der Geräterückseite ( $\rightarrow$  Abb. 5, Seite 21) über das mitgelieferte Netzkabel mit einer geerdeten Steckdose. Eine manuelle Anpassung an die länderspezifische Netzspannung ist nicht erforderlich.

**Warnung:** Verwenden Sie nur die für das Gerät bereitgestellten Netzkabel.

Verwenden Sie keine Mehrfachsteckdosen oder Verlängerungskabel. Die Verwendung von defekten Mehrfachsteckdosen oder Verlängerungskabeln kann zu Personenschäden oder Schäden am Gerät führen.

### 3.4.4 Anschluss des Signalkabels (Digital I/O)

Verwenden Sie zum Anschluss von externen Geräten an die Digital I/O-Buchsen auf der Geräterückseite das 6-adrige Mini-DIN-Signalkabel (Best.-Nr. 6000.1004).

- Verbinden Sie den 6-poligen Stecker des Kabels mit dem Port Digital I/O 1 (2, 3 oder 4). Informationen zur Anschlussbelegung der Buchsen und des Kabels finden Sie auf der Seite 191.
- Verbinden Sie f
  ür jeden Relaisausgang und digitalen Eingang, den Sie verwenden möchten, die Signal- und Erdungsadern des Kabels mit den entsprechenden Anschl
  üssen des externen Ger
  ätes (→ Dokumentation f
  ür das externe Ger
  ät).

**Hinweis:** Folgen Sie zum Anschluss einer Pumpe des Typs LPG-3400XRS den Schritten in der *Bedienungsanleitung* zur Pumpe.

- 3. Nehmen Sie bei der Konfiguration des Autosamplers im Programm **Server Configuration** von Chromeleon folgende Einstellungen vor:
  - Wählen Sie auf der Seite **Relays** den entsprechenden Relaisausgang aus (→ Seite 44).
  - Wählen Sie auf der Seite Inputs den entsprechenden digital Eingang aus (→ Seite 44).

## 3.5 Einrichten des Autosamplers in Chromeleon

Die nachfolgenden Seiten geben Ihnen einen kurzen Überblick, wie Sie den Autosampler in Chromeleon einrichten. Einzelheiten hierzu finden Sie in der *Chromeleon-Hilfe*. Informationen zum Einrichten des Autosamplers unter DCMSLink finden Sie in Kapitel 3.6 ( $\rightarrow$  Seite 46).

**I** Hinweis: Wenn der Autosampler mit dem Chromeleon-Rechner verbunden ist, vergewissern Sie sich, dass die Chromeleon-Software auf dem Rechner installiert ist, *bevor* Sie den Autosampler zum ersten Mal einschalten. Nur dann wird der USB-Treiber für den Autosampler automatisch geladen und das Windows<sup>®</sup>-Betriebssystem erkennt den Autosampler automatisch, wenn dieser eingeschaltet wird.

#### 3.5.1 Laden des USB-Treibers

- 1. Schalten Sie, sofern noch nicht geschehen, den Chromeleon-Rechner ein.
- Melden Sie sich unter Windows Vista<sup>®</sup> (Windows<sup>®</sup> XP, Windows<sup>®</sup> 7 oder Windows<sup>®</sup> Server 2008) als
  - Administrator an, wenn es sich um einen lokalen Rechner handelt.
  - Anwender mit Administratorrechten an, wenn es sich um einen Netzwerkrechner handelt.
- Öffnen Sie das Programm Chromeleon Server Monitor über einen Doppelklick auf das Server Monitor-Symbol in der Windows Taskleiste.

Wenn das Server Monitor-Symbol nicht in der Taskleiste vorhanden ist, klicken Sie auf **Start** in der Taskleiste. Wählen Sie **Programme** (oder **Alle Programme**, abhängig vom Betriebssystem), wählen Sie dann **Chromeleon**, und klicken Sie danach auf **Server Monitor**.

- 4. Klicken Sie auf **Start**, um den Server zu starten.
- Schließen Sie das Server Monitor-Fenster mit Close. Das Server Monitor-Symbol serscheint in der Taskleiste.
  - **Hinweis:** Über die Schaltfläche **Quit Monitor** können Sie das Programm **Server Monitor** verlassen, der Server wird jedoch nicht gestoppt. Um den Server anzuhalten, klicken Sie auf **Stop**.
- 6. Schalten Sie den Autosampler über den Netzschalter auf der Rückseite ein.
- 7. *Windows Vista, Windows 7 und Windows Server 2008* erkennen den neuen Autosampler automatisch und führen die USB-Installation durch.

Wird der Autosampler nicht automatisch erkannt und stattdessen ein Installationsassistent gestartet, deutet dies darauf hin, dass Sie den Autosampler mit dem Rechner verbunden und eingeschaltet haben, ohne dass die Chromeleon-Software installiert ist.

Gehen Sie folgendermaßen vor:

- a) Brechen Sie den Assistenten ab.
- b) Schalten Sie den Autosampler aus.
- c) Installieren Sie Chromeleon.
- d) Schalten Sie den Autosampler ein. Windows erkennt jetzt den Autosampler und installiert die USB-Software für den Autosampler automatisch.
- 8. Windows XP

erkennt den neuen Autosampler automatisch und startet einen Assistenten (**Neue Hardware gefunden**), der Sie durch die weitere USB-Installation führt. Wählen Sie die folgenden Optionen:

- a) Falls eine Meldung erscheint, ob eine Verbindung mit Windows Update hergestellt werden soll, um nach Software zu suchen, wählen Sie **Nein, diesmal nicht**.
- b) Akzeptieren Sie die standardmäßig ausgewählte Option **Software automatisch installieren** und klicken Sie auf **Weiter**>.
- c) Klicken Sie auf **Fertigstellen**, wenn der Assistent meldet, dass die Software für den Autosampler installiert wurde.

Wenn Windows den Autosampler nicht automatisch erkennt und stattdessen nach einer USB-Konfigurationsdatei (cmwdmusb.inf) fragt, deutet dies darauf hin, dass Sie den Autosampler mit dem Rechner verbunden und eingeschaltet haben, ohne dass die Chromeleon-Software installiert ist. Gehen Sie folgendermaßen vor:

- a) Klicken Sie in der Meldung von Windows auf Abbrechen.
- b) Schalten Sie den Autosampler aus.
- c) Installieren Sie Chromeleon.
- d) Schalten Sie den Autosampler ein. Windows erkennt den Autosampler jetzt automatisch und startet den Assistenten **Neue Hardware** gefunden.

#### 3.5.2 Installieren des Autosamplers

Nachdem die USB-Software für den Autosampler installiert ist ( $\rightarrow$  Seite 37), können Sie den Autosampler in Chromeleon installieren und konfigurieren:

- 1. Starten Sie, falls erforderlich, den Chromeleon Server Monitor und den Chromeleon Server ( $\rightarrow$  Seite 37).
- Starten Sie das Installationsprogramm Server Configuration von Chromeleon. Klicken Sie auf Start in der Taskleiste. Wählen Sie Programme (oder Alle Programme, abhängig vom Betriebssystem), wählen Sie dann Chromeleon und klicken Sie danach auf Server Configuration.
- 3. Klicken Sie, falls erforderlich, auf das Pluszeichen neben dem Servernamen, um sich die Einträge unterhalb des Servers anzeigen zu lassen.
- 4. Wählen Sie die Zeitbasis aus, der Sie den Autosampler zuordnen möchten, oder legen Sie eine neue Zeitbasis an (über Add Timebase im Menü Edit).
- 5. Öffnen Sie das Dialogfenster Add device to timebase. Wählen Sie dazu im Menü Edit den Befehl Add Device (oder wählen Sie die Zeitbasis per Rechtsklick aus und wählen Sie im Menü den Punkt Add Device).
- 6. Wählen Sie unter Manufacturers den Eintrag Dionex HPLC: UltiMate 3000 und dann unter Devices den Eintrag WPS-3000(RS) Autosampler aus.
- 7. Ein Assistent führt Sie durch die Installation. Chromeleon stellt eine Verbindung zum Autosampler her und übernimmt die Einstellungen der Gerätefirmware nach Chromeleon. Bestätigen Sie die Meldung mit OK.
- Prüfen Sie die Einstellungen auf den einzelnen Seiten auf Richtigkeit und nehmen Sie gegebenenfalls weitere Einstellungen vor. Die einzelnen Seiten sind im Kapitel 3.5.3.1 beschrieben (→ Seite 40).
- 9. Klicken Sie auf **Fertig stellen**, um die Basiskonfiguration des Autosamplers zu beenden.
- 10. Speichern Sie die Konfiguration mit **Save Installation** im Menü **File** und schließen Sie das Serverkonfigurationsprogramm.

#### 3.5.3 Konfigurieren des Autosamplers

#### 3.5.3.1 Erstinstallation

Bei der Installation des Autosamplers werden die Einstellungen der Gerätefirmware nach Chromeleon übertragen. Prüfen Sie diese Standardeinstellungen auf Richtigkeit und nehmen Sie gegebenenfalls weitere Einstellungen vor. Sie können die Konfigurationsseiten auch nach der Installation erneut öffnen, wenn Sie die Einstellungen verändern möchten ( $\rightarrow$  Seite 45).

**Hinweis:** Änderungen, die Sie im Dialogfenster **Commands**, in einem Programm (PGM) oder auf einem Steuerfenster vornehmen, haben keinen Einfluss auf die Standardeinstellungen auf diesen Registerkarten.

Weitere Informationen zu einer Seite erhalten Sie über die Schaltfläche Hilfe.

#### Seite General

Die Seite General zeigt die allgemeinen Geräteparameter:

General		×
Device Name:	Sampler	
Demo <u>M</u> ode	Firmware Down <u>l</u> oad	
Module <u>A</u> ddress:		
		Browse

Abb. 9: Seite General

#### • Device Name

zeigt den Namen an, unter dem der Autosampler in der Installationsumgebung und im Chromeleon Client geführt wird. Ändern Sie den vorgegebenen Namen nicht, wenn Sie den Autosampler über vorhandene Steuerfenster steuern möchten. Wenn Sie einen anderen Namen eingeben, müssen Sie gegebenenfalls die Links der Bedienelemente auf den Steuerfenstern und den Namen des Autosamplers in den Programmen entsprechend anpassen.

#### • Demo Mode

Achten Sie darauf, dass dieses Kontrollkästchen nicht ausgewählt ist. Im Demo Mode simuliert Chromeleon die Funktionen des Autosamplers. Wenn der Demo Mode eingeschaltet ist, kann im Feld **Module Address** keine Moduladresse ausgewählt werden. Wenn Sie die Seite **General** verlassen, ohne eine Moduladresse eingegeben zu haben, wird automatisch der Demomodus aktiviert.

#### • Module Address

Wählen Sie die Moduladresse des Autosamplers aus. Klicken Sie dazu auf die Schaltfläche **Browse** und wählen Sie den Autosampler per Doppelklick aus. Seine Adresse wird damit automatisch in das Feld **Module Address** übernommen.

#### • Firmware Download

Über diese Schaltfläche kann die Firmware-Version, die in Chromeleon für den Autosampler enthalten ist, von Chromeleon auf den Autosampler übertragen werden. (Die Schaltfläche ist ausgeblendet, wenn der Demo Mode eingeschaltet ist.)

Die aktuelle Firmware-Version ist bei Auslieferung des Autosamplers installiert. Sollte dennoch eine Aktualisierung der Firmware erforderlich sein, folgen Sie der Beschreibung im Kapitel 7.10 ( $\rightarrow$  Seite 159).

Sobald Sie bei der Erstinstallation die Seite **General** (bei ausgeschaltetem Demo Mode) verlassen, versucht Chromeleon, eine Verbindung zum Autosampler herzustellen und die Einstellungen der Gerätefirmware nach Chromeleon zu übernehmen und nimmt dann die entsprechenden Einstellungen auf den folgenden Seiten des Assistenten vor. Bestätigen Sie die entsprechende Meldung mit **OK**.

#### Seite Sharing

Standardmäßig ist der Autosampler der Zeitbasis zugeordnet, unter der er installiert wird. Sie können den Autosampler jedoch auch auf zwei Zeitbasen gemeinsam nutzen. Wählen Sie dazu eine zweite Zeitbasis aus der Liste der verfügbaren Zeitbasen aus. Der Autosampler erscheint dann in der Serverkonfiguration unter den beiden Zeitbasen und kann von beiden Zeitbasen aus gesteuert werden.

#### Seite Segments / Pump Link

Auf dieser Seite wird angezeigt, welche Probengefäße standardmäßig in welchem Segment installiert sind. Wenn Ihr UltiMate 3000-System eine Pumpe aus der UltiMate 3000-Pumpenserie enthält, sollten Sie hier außerdem den Autosampler mit der Pumpe verbinden. Nehmen Sie diese Einstellung grundsätzlich vor, da sie:

- angibt, welche Pumpe den Fluss liefert. Diese Information ist wichtig, wenn Sie im Bypass-Modus arbeiten (→ Seite 13), da Chromeleon bei der Berechnung, wann das Injektionsventil in die Position Load umgeschaltet wird, die Flussrate berücksichtigt.
- erlaubt, den Injektionsbefehl des Autosamplers mit den Kolbenhüben einer Niederdruckgradientenpumpe DGP-3600 bzw. LPG-3400 zu synchronisieren. Damit wird sichergestellt, dass alle Injektionen zur selben Phase des Pumpenzyklus erfolgen. Auf diese Weise werden die analytischen Ergebnisse deutlich verbessert.

Segments / Pump Link		
Wel	Plate Sampler Segment Types	
Re	d 40 Standard Vials	
Gn	een 40 Standard Vials 💌	
Blu	ue 40 Standard Vials 💌	
W	aming: Ensure that these settings match the installed items!	
- Pum Flo	p Link w through <u>S</u> ampler is delivered by pump (s):	
<none></none>		
<none></none>		
UI	M3PUMP_L_STRK M3PUMP_R_STRK pump is an EPG-3x00 or a DGP-3600.	

Abb. 10: Seite Segments / Pump Link (hier: wenn das UltiMate 3000-System eine Pumpe DGP-3600 enthält)

#### • Well Plate Sampler Segment Types

Zeigt an, welche Probengefäße standardmäßig im roten, grünen und blauen Segment installiert sind. Prüfen und ändern Sie diese Einstellungen, falls erforderlich. Wenn Sie zu einem späteren Zeitpunkt andere Probengefäße im Autosampler installieren, müssen Sie die Einstellung in Chromeleon entsprechend anpassen ( $\rightarrow$  Seite 84).

#### • Pump Link

Geben Sie an, mit welcher Pumpe eines UltiMate 3000-Systems der Autosampler standardmäßig verbunden werden soll. Wählen Sie die Pumpe aus der Liste **Flow through sampler is delivered by pump(s)** aus. Die Liste zeigt alle Pumpen, die auf der Zeitbasis installiert sind.

**Hinweis:** Wenn es sich bei der Pumpe um eine LPG-3400XRS handelt *oder* wenn Sie eine LPG-3400XRS *und* eine weitere Pumpe mit dem Autosampler verbinden möchten, folgen Sie bitte der Beschreibung in der *Bedienungsanleitung zur LPG-3400XRS*.

- Wenn das UltiMate 3000-System eine Pumpe des Typs DGP-3600 enthält Wählen Sie UM3PUMP\_L\_STRK, wenn der Autosampler standardmäßig mit der linken Pumpe verbunden werden soll. Wählen Sie UM3PUMP\_R\_STRK, wenn der Autosampler standardmäßig mit der rechten Pumpe verbunden werden soll.
- *Wenn das UltiMate 3000-System eine andere Pumpe als eine DGP-3600 enthält* Wählen Sie UM3PUMP\_STROKE.
- Wenn der Autosampler nicht mit einer Pumpe verbunden werden soll Wählen Sie <None>.

Nur wenn Sie auf dieser Seite den Autosampler mit einer DGP-3600 oder LPG-3400 verbunden haben, ist die Synchronisierung des Injektionsbefehls mit den Kolbenhüben der Pumpe möglich.

Im Chromeleon Client können Sie dann für die jeweilige Anwendung festlegen, ob Sie mit oder ohne Synchronisierung arbeiten möchten und gegebenenfalls eine andere Pumpe festlegen ( $\rightarrow$  Seite 82).

#### **Seite Options**

Auf dieser Seite wird angezeigt, für welche Anwendung der Autosampler konfiguriert ist oder ob eine anwenderspezifische Konfiguration vorhanden ist. Prüfen und ändern Sie gegebenenfalls die angezeigte Konfiguration.

O <u>M</u> icro	C <u>C</u> ustom	
Analytical	Sample Joop volume (µl):	130
O Semiprep	Syringe size (µl):	100
C Low Dispersion	Seat capillary <u>v</u> olume (µl):	3.1
✓ Thermostatting option installed		

Abb. 11: Seite Options

#### • Micro, Analytical, Semiprep, Low Dispersion und Custom

Micro, Analytical, Semiprep und Low Dispersion zeigen an, für welche Anwendung der Autosampler vorkonfiguriert ist. Die genaue Konfiguration wird unterhalb von Custom angezeigt. Diese Angaben dienen der Information und können nicht verändert werden. Klicken Sie Custom an, wenn Sie stattdessen eine spezifische Konfiguration verwenden möchten, und machen Sie die entsprechenden Angaben in den Feldern unterhalb von Custom. Beachten Sie bitte auch die Informationen für die Einrichtung eines WPS-3000TBRS und WPS-3000TBSL Analytical weiter unten.

#### Sample loop volume (μL)

Volumen der Probenschleife (= Gradientenverzögerungsvolumen) Geben Sie bei Dionex-Probenschleifen den Wert ein, der auf dem Aufkleber der Probenschleife in Klammern vermerkt ist (V =  $xx \mu L$ ).

- Syringe size (μL) Spritzengröße
- Seat capillary volume (μL) Volumen der Nadelsitzkapillare

#### WPS-3000TBRS

Wenn Sie einen WPS-3000TBRS einrichten, ist **Custom** ausgewählt. Für die Standardkonfiguration des Autosamplers sind keine Änderungen erforderlich. Wenn Sie jedoch anstelle der 25- $\mu$ L-Probenschleife die 100- $\mu$ L-Probenschleife aus dem Zubehör installieren, müssen Sie den Wert für Sample **loop volume** auf **130** ändern.

#### WPS-3000TBSL Analytical

Wenn Sie einen WPS-3000TBSL Analytical einrichten möchten, müssen Sie den Wert für **Sample loop volume** anpassen. Klicken Sie **Custom** an und geben Sie für **Sample loop volume** 150 ein.

#### • Thermostatting option installed

Wenn es sich bei Ihrem Autosampler um ein Gerät mit Probenthermostatisierung handelt, vergewissern Sie sich, dass das Kontrollkästchen angeklickt ist.

#### • Automatically Wash after Abort errors

Das Kontrollkästchen ist standardmäßig aktiviert. Übernehmen Sie diese Einstellung, wenn Chromeleon nach einem Abortfehler automatisch einen Waschzyklus durchführen soll.

Wenn der Autosampler mit Chromeleon verbunden wird ("Connect"), werden die Einstellungen auf den Seiten **Segments / Pump Link** und **Options** mit der Gerätekonfiguration verglichen, die der Autosampler beim Aufbau der Verbindung an Chromeleon meldet. Stimmen Gerätekonfiguration und Einstellungen in Chromeleon nicht überein, wird der Autosampler von Chromeleon getrennt ("Disconnect") und es erscheint eine Meldung im Chromeleon Audit Trail. Prüfen Sie die Einstellungen und gleichen Sie diese gegebenenfalls über die Schaltflächen **Retrieve configuration from module** oder **Send configuration to module** auf der Registerkarte **General** ab ( $\rightarrow$  Seite 45).

#### Seite Relays

Die Seite **Relays** listet alle zur Verfügung stehenden Relais auf (Relay\_1 bis Relay\_4). Klicken Sie ein Kontrollkästchen an, um das entsprechende Relais zu aktivieren. Nicht aktivierte Relais stehen unter Chromeleon nicht zur Verfügung. Sie können den Namen und die Zeitbasis-Zuordnung für ein Relais ändern. Klicken Sie dazu den Relaisnamen doppelt an (oder drücken Sie auf der Tastatur **F2**) und nehmen Sie die Änderungen in der Dialogbox **Device Configuration** vor.

#### **Seite Inputs**

Die Seite **Inputs** listet alle zur Verfügung stehenden Digitaleingänge auf (Input\_1 bis Input\_4). Klicken Sie ein Kontrollkästchen an, um den entsprechenden Digitaleingang zu aktivieren. Nicht aktivierte Eingänge stehen unter Chromeleon nicht zur Verfügung. Sie können den Namen und die Zeitbasis-Zuordnung für einen Digitaleingang ändern. Klicken Sie dazu den Name des Eingangs doppelt an (oder drücken Sie auf der Tastatur **F2**) und nehmen Sie die Änderungen in der Dialogbox **Device Configuration** vor.

#### 3.5.3.2 Ändern der Konfiguration

Sie können die Konfigurationsseiten auch später erneut öffnen, wenn Sie die Standardeinstellungen verändern möchten.

- 1. Starten Sie das Programm Server Configuration ( $\rightarrow$  Seite 39).
- 2. Markieren Sie den Autosampler in der Zeitbasis mit einem Rechtsklick und wählen Sie im Menü den Punkt **Properties**.
- 3. Ändern Sie die Einstellungen auf den Registerkarten wie gewünscht ab. Die einzelnen Seiten sind im Kapitel 3.5.3.1 beschrieben (→ Seite 40).

Auf der Registerkarte **General** stehen nun zwei weitere Schaltflächen zur Verfügung, mit denen Sie die Konfiguration des Autosamplers und die Einstellungen in Chromeleon abgleichen können:

Schaltfläche	Beschreibung
Retrieve configuration from module	Drücken Sie diese Schaltfläche, um die aktuelle Konfiguration des Autosamplers nach Chromeleon zu übertragen. Die Einstellungen auf den Registerkarten <b>Segments / Pump Link</b> und <b>Options</b> werden in Chromeleon aktualisiert. Dies ist jedoch nur erforderlich, wenn Sie einen anderen Autosampler anschließen oder wenn die Automatik beim Start des Assistenten fehlschlägt.
Send configuration to module	Drücken Sie diese Schaltfläche, um die aktuell in Chromeleon auf den Registerkarten <b>Segments / Pump Link</b> und <b>Options</b> festgelegte Konfiguration zum Autosampler zu übertragen.

Zusätzlich zu den Seiten des Installationsassistenten steht die Registerkarte **Error Levels** zur Verfügung. Auf dieser Registerkarte ist festgelegt, welchen Stellenwert die einzelnen Meldungen unter Chromeleon haben. In der Regel sind hier keine Änderungen erforderlich.

4. Speichern Sie die geänderte Konfiguration mit **Save Installation** im Menü **File** und schließen Sie das Serverkonfigurationsprogramm.

## 3.6 Einrichten des Autosamplers in DCMSLink

- 1. Installieren und konfigurieren Sie die DCMSLink-Software (→ *DCMSLink Installation Guide*). Der Guide steht auf der DCMSLink-DVD im Verzeichnis *Additional Documents*\*DCMSLink User Documents* zur Verfügung.
- 2. Öffnen Sie das Programm **Server Configuration** von Chromeleon ( $\rightarrow$  *DCMSLink Installation Guide*).
- 3. Nehmen Sie den Autosampler im Programm Server Configuration in eine Zeitbasis auf. Die Vorgehensweise entspricht der Beschreibung in Kapitel 3.5.2 (→ Seite 39).
- 4. Konfigurieren Sie den Autosampler. Die Vorgehensweise entspricht der Beschreibung in Kapitel 3.5.3 (→ Seite 40).

Weitere Informationen zu DCMSLink finden Sie in der *DCMSLink-Hilfe* und im *DCMSLink Quick Start Guide*. Der Quick Start Guide steht ebenfalls auf der DCMSLink-DVD zur Verfügung.

## 4 Vorbereitung für den Betrieb (Inbetriebnahme)

## 4.1 Übersicht

# ▲ Vorsicht: Der Autosampler ist bei Auslieferung mit einer 2-Propanol-Lösung gefüllt. Achten Sie bei der ersten Inbetriebnahme darauf, dass das verwendete Lösungsmittel mit 2-Propanol mischbar ist. Ist dies nicht der Fall, stellen Sie das Lösungsmittel schrittweise um.

Achten Sie beim Anschluss von Kapillaren darauf, dass keine Verunreinigungen an den Anschlüssen haften. Eindringende Schmutzpartikel können zu Schäden am System führen.

Nachdem Sie den Autosampler ausgepackt, aufgestellt und angeschlossen haben ( $\rightarrow$  Kapitel 3.1 bis 3.4, ab Seite 31), bereiten Sie den Autosampler für den Betrieb vor:

- 1. Schließen Sie den Autosampler fluidisch an die Pumpe bzw. den Säulenthermostaten oder die Trennsäule an.
- 2. Prüfen Sie die Anschlüsse am Injektionsventil ( $\rightarrow$  Seite 27).
- 3. Schließen Sie die Waschflüssigkeit an ( $\rightarrow$  Seite 50).
- 4. Prüfen Sie die Anschlüsse am Spritzenventil ( $\rightarrow$  Seite 53).
- 5. Schließen Sie die Drainage an ( $\rightarrow$  Seite 54).
- 6. Wenn Sie den Autosampler mit Chromeleon betreiben möchten Installieren Sie den Autosampler in Chromeleon ( $\rightarrow$  Seite 37).
- 7. Schalten Sie den Autosampler ein ( $\rightarrow$  Seite 63).
- 8. Spülen Sie die Spritze ( $\rightarrow$  Seite 55).
- 9. Installieren und bestücken Sie die Probenhalter oder Mikrotiterplatten ( $\rightarrow$  Seite 57).
- 10. Bestücken Sie die 5-Positionen Halter ( $\rightarrow$  Seite 59).
- Nur Autosampler mit Probenthermostatisierung Schalten Sie die Probenthermostatisierung ein und legen Sie die Temperatur fest (→ Seite 83).
- Prüfen und ändern Sie gegebenenfalls die Einstellung für die Leakerkennung (→ Seite 86).

- Passen Sie, falls erforderlich, die Helligkeit und den Kontrast der Displayanzeige an Ihre Anforderungen an (→ Seite 86).
- 14. Ehe Sie mit der Probenanalyse beginnen, äquilibireren Sie das gesamte System  $(\rightarrow$  Seite 60).

## 4.2 Hinweise zum Anschluss von Kapillaren

Beachten Sie beim Anschluss von Kapillaren die folgenden Hinweise:

- Beachten Sie die Sicherheitshinweise zu Kapillaren und Kapillarverbindungen im Kapitel 1.2.2 (→ Seite 4).
- Achten Sie darauf, dass keine Verunreinigungen an den Anschlüssen haften. Eindringende Schmutzpartikel können zu Schäden am System führen.
- Innerhalb eines UltiMate 3000-Systems werden unterschiedliche Fittingsysteme verwendet. Installieren Sie die Kapillarverbindungen nur an der dafür vorgesehenen Position.
- Verwenden Sie nur die mitgelieferten, fertig konfektionierten Anschlusskapillaren bzw. Original-Dionex-Ersatzkapillaren.
- Verwenden Sie möglichst Viper-Kapillarverbindungen. Beachten Sie die Hinweise in der Anleitung, die der Kapillare beiliegt.
- Abhängig von der verwendeten Fittingverbindung, beachten Sie außerdem die folgenden Punkte:
  - Viper-Fittingverbindungen

Lösen oder ziehen Sie Viper-Fittingverbindungen *nur* mit der Rändelschraube und *nur* per Hand fest (verwenden Sie *kein* Werkzeug).

Ziehen Sie die Schraube zunächst handfest an. Wenn an der Verbindung eine Undichtigkeit auftreten sollte, ziehen Sie die Schraube etwas weiter fest. Bleibt die Undichtigkeit bestehen, entfernen Sie die Kapillare, reinigen Sie die Kapillarenden vorsichtig mit einem mit Isopropanol getränkten Tuch und bauen Sie die Kapillare wieder ein. Verwenden Sie eine andere Viper-Kapillare, wenn die Undichtigkeit weiterhin bestehen bleibt.

**Hinweise:** Kapillaren mit Viper-Fittingen können Sie für unterschiedliche Verbindungen wieder verwenden.

Schwarze Rändelschrauben (wie in Abb. 12 gezeigt) können Sie von der Kapillare entfernen und wieder aufsetzen.



Abb. 12: Viper-Fittingverbindung (Beispiel)

#### • Herkömmliche (nicht-Viper) Fittingverbindungen

Ziehen Sie diese Fittingverbindungen nicht zu fest an. Ziehen Sie die Verbindung gegebenenfalls nach, wenn eine Undichtigkeit auftritt.

Bleibt die Undichtigkeit bestehen, reinigen Sie zunächst den Anschlussport mit einem Reinigungsstäbchen (Best.-Nr. 6040.0006) säubern. Wechseln Sie die Kapillare und/oder das Fitting, wenn die Undichtigkeit weiterhin bestehen bleibt. Um ein erhöhtes Totvolumen oder Beschädigungen und Undichtigkeiten zu vermeiden, verwenden Sie bereits benutzte Fittingverbindungen nur für dieselbe Kapillarverbindung wieder.

## 4.3 Anschließen der Waschflüssigkeit

Thermo Fisher Scientific empfiehlt, die Waschflüssigkeit kontinuierlich zu entgasen. Wenn das UltiMate 3000-System eine Pumpe LPG-3400 oder ein geeignetes SRD-3x00 Solvent Rack enthält, können Sie die Waschflüssigkeit online über den Degaser der Pumpe oder des Solvent Racks entgasen ( $\rightarrow$  Kapitel 4.3.1).

Wenn dies nicht möglich ist, z.B. weil Ihr UltiMate 3000-System eine semipräparative Pumpe HPG-3200P enthält, können Sie alternativ die Waschflasche aus dem Zubehör installieren  $(\rightarrow$  Seite 52).

#### 4.3.1 Anschluss der Waschflüssigkeit bei Online-Entgasung

Schließen Sie die Waschflüssigkeit wie folgt an den Degaser an:

- Direkt, wenn ein freier Degaserkanal zur Verfügung steht (siehe unten) In diesem Fall können Sie die Waschflüssigkeit frei wählen.
- Indirekt, über die Pumpe des UltiMate 3000 Systems ( $\rightarrow$  Seite 51) Bei dieser Methode muss eines der Laufmittel der Pumpe als Waschlösung geeignet sein.

Alle für den Anschluss erforderlichen Teile sind im Online Degas Wash Kit enthalten. Das Kit ist Teil des Zubehörs.

#### Direkter Anschluss an den Degaser

- 1. Füllen Sie eine Vorratsflasche mit einer geeigneten Waschflüssigkeit.
- Bereiten Sie zunächst den Ansaugschlauch für die Waschflüssigkeit vor und platzieren 2. Sie ihn in die Vorratsflasche. Verbinden Sie dann die Vorratsflasche mit dem Degasereinlass. (Die Vorgehensweise entspricht dem Anschluss des Eluentenvorrats. Einzelheiten finden Sie unter Anschluss des Eluentenvorrats in der Bedienungsanleitung zur Pumpe bzw. zum Solvent Rack des UltiMate 3000-Systems.)
- 3. Stellen Sie die Vorratsflasche in die Wanne des Solvent Racks.
- 4 Schließen Sie den langen Schlauch aus dem Online Degas Wash Kit am Anschlussport für die Waschflüssigkeit am Spritzenventil an.



Abb. 13: Anschluss der Waschflüssigkeit am Spritzenventil

5. Kürzen Sie das freie Schlauchende mit einem geeigneten Kapillarschneider entsprechend Ihren Anforderungen. Achten Sie auf einen rechtwinkligen Schnitt.

- 6. Verbinden Sie das freie Schlauchende mit dem Degaserauslass. Verwenden Sie dazu die Rändelschraube und das Fitting aus dem Online Degas Wash Kit.
- 7. Entlüften Sie die Spritze, bis keine Luftblasen mehr sichtbar sind ( $\rightarrow$  Seite 55).

#### Indirekter Anschluss an den Degaser über die Pumpe des UltiMate 3000-Systems

- 1. Füllen Sie eine Vorratsflasche mit einer geeigneten Waschflüssigkeit.
- Bereiten Sie zunächst den Ansaugschlauch für die Waschflüssigkeit vor und platzieren Sie ihn in der Vorratsflasche. Verbinden Sie dann die Vorratsflasche mit dem Degasereinlass. (Die Vorgehensweise entspricht dem Anschluss des Eluentenvorrats. Einzelheiten finden Sie unter Anschluss des Eluentenvorrats in der Bedienungsanleitung zur Pumpe bzw. zum Solvent Rack des UltiMate 3000-Systems.)
- 3. Stellen Sie die Vorratsflasche in die Wanne des Solvent Racks.
- 4. Schließen Sie den langen Schlauch aus dem Online Degas Wash Kit am Anschlussport für die Waschflüssigkeit am Spritzenventil an (→ Abb. 13, Seite 50).
- 5. Verbinden Sie das andere Ende dieses Schlauchs mit dem T-Stück aus dem Online Degas Wash Kit.
- 6. Lösen Sie die Verbindung zwischen dem Degaserauslass und der Pumpe, je nach Pumpentyp am Sternventil oder am Pumpeneinlass. Schließen Sie den Schlauch, der vom Degaserauslass kommt, an das T-Stück an.
- 7. Stellen Sie die Verbindung zwischen dem T-Stück und dem Sternventil der Pumpe bzw. dem Pumpeneinlass her. Verwenden Sie dazu den kurzen Schlauch aus dem Online Degas Wash Kit.



Anschluss des Schlauchs zur Pumpe (hier zum Sternventil)

Anschluss des Schlauchs vom Degaserauslass

T-Stück

Abb. 14: Beispiel für den Anschluss der Waschflüssigkeit über die Pumpe

8. Entlüften Sie die Spritze, bis keine Luftblasen mehr sichtbar sind ( $\rightarrow$  Seite 55).

Anschluss des Schlauchs

vom Spritzenventil

#### 4.3.2 Anschluss der Waschflasche

Wenn eine Online-Entgasung der Waschflüssigkeit innerhalb des UltiMate 3000-Systems nicht möglich ist, können Sie alternativ die 125-mL-Waschflasche aus dem Zubehör verwenden. Die Waschflasche wird im Probenraum des Autosamplers installiert. Der Halteclip der Waschflasche ( $\rightarrow$  Abb. 16) ist bereits vorinstalliert.

- 1. Füllen Sie die Flasche mit einer geeigneten Waschflüssigkeit. Entgasen Sie die Waschflüssigkeit vor Verwendung im Ultraschallbad mit Unterdruck oder entgasen Sie die Flüssigkeit kontinuierlich mit einem Vakuumfiltrationsgerät.
- 2. Prüfen Sie, ob der Schlauch für die Waschflüssigkeit durch die Öffnung im Deckel ausreichend tief in die Waschflasche geführt ist, möglichst bis zum Boden. (Halten Sie die Flasche gegebenenfalls gegen das Licht.)
- 3. Schließen Sie das andere Ende des Schlauchs am Spritzenventil an ( $\rightarrow$  Abb. 15).
  - **Hinweis:** Um zu verhindern, dass Luft in das System gelangt, muss der Schlauch beim Anschließen an das Spritzenventil komplett mit der Waschflüssigkeit gefüllt sein. Drücken Sie die Waschflasche zusammen und verschließen Sie dabei das Belüftungsloch im Flaschendeckel mit dem Daumen. So wird Druck in der Flasche aufgebaut und der Schlauch wird mit Flüssigkeit gefüllt.



Anschlussport Waschflasche

Abb. 15: Anschluss der Waschflasche am Spritzenventil

4. Hängen Sie die Halteklammer mit der Waschflasche in den Halter ein. Die Halteklammer muss im Halter einrasten.







Waschflasche mit Halteklammer (Rückansicht)

Halter Waschflasche

Abb. 16: Einsetzen der Waschflasche

5. Entlüften Sie Spritze, bis keine Luftblasen mehr sichtbar sind ( $\rightarrow$  Seite 55).

Um die Waschflasche neu zu füllen oder die Waschflüssigkeit auszutauschen, greifen Sie die Flasche einschließlich der Halteklammer und schieben Sie beides senkrecht nach oben (nicht verkanten). Die Halteklammer rastet aus. Ziehen Sie die Flasche einschließlich der Halteklammer heraus.

**Hinweis:** Wenn Sie für Ihre Applikation mehr Waschflüssigkeit benötigen, können Sie einen längeren Schlauch (mit Flansch für das Ventilfitting) anschließen und eine größere Waschflasche neben den Autosampler oder in die Wanne des Solvent Racks stellen. (Sie können beispielsweise den Schlauch aus dem Online Degas Wash Kit verwenden, das im Zubehör enthalten ist.) Um diesen längeren Schlauch vollständig zu füllen, müssen Sie den Spülvorgang unter Umständen mehrfach wiederholen.

## 4.4 Anschlüsse am Spritzenventil

Prüfen Sie am Spritzenventil, ob die Pufferschleife und die Spritze sowie die Schläuche für die Versorgung des Waschports und die Waschflüssigkeit korrekt angeschlossen sind. Ziehen Sie die Verbindungen gegebenenfalls nach.



Abb. 17: Anschlüsse am Spritzenventil

Wenn Sie die Verbindung der Spritze oder Waschflüssigkeit nachgezogen haben ( $\rightarrow$  Seite 55), spülen Sie die Spritze, um Luftblasen aus der Spritze, dem Spritzenventil und den Ansaugschläuchen für die Reinigungslösung zu entfernen.

Wenn Sie die Verbindung der Pufferschleife nachgezogen haben, führen Sie einen Waschzyklus für die Pufferschleife durch, um Luftblasen aus der Pufferschleife zu entfernen. Führen Sie in Chromeleon den Befehl **WashBufferLoop** aus oder wählen Sie am Gerätedisplay im Menü **Control** den Befehl **Buffer loop wash**.

## 4.5 Anschließen der Drainage

Zur Ableitung von Flüssigkeiten aus dem Geräteinneren verfügt der Autosampler rechts unterhalb des Gerätes über *zwei* oder *drei* Abläufe (je nach Version).

- Der Autosampler ohne Probenthermostatisierung verfügt über *zwei* Abläufe, über die der Spritzen- und Waschabfall sowie eventuelle Leckage-Flüssigkeiten abgeleitet werden.
- Der Autosampler mit Probenthermostatisierung verfügt über *drei* Abläufe. Über die vorderen Abläufe werden der Spritzen- und Waschabfall sowie eventuelle Leckage-Flüssigkeiten abgeleitet. Der hintere Ablauf dient zur Ableitung von Kondenswasser aus dem Drehteller.



Abb. 18: Drainage-Abläufe

Leiten Sie die Flüssigkeiten über das Drainagesystem des UltiMate 3000-Systems in den Abfall. Die entsprechenden Komponenten stehen im Drainage-Kit für das UltiMate 3000 System zur Verfügung. Das Kit ist im Lieferumfang der UltiMate 3000-Pumpen enthalten und kann auch separat bestellt werden (Best.-Nr. 6040.0005). Es enthält alle erforderlichen Komponenten für die Systemdrainage sowie eine detaillierte Installationsanleitung.

## 4.6 Entfernen von Luftblasen aus der Spritze (Prime)

Luft in der Spritze kann die Messergebnisse stark beeinflussen. Wählen Sie eine der unten genannten Alternativen, um die Spritze zu spülen und die Luft aus der Spritze zu entfernen.

- Spülen Sie die Spritze über Chromeleon.
- Spülen Sie die Spritze vom Gerätedisplay aus.

Alle mit dem Spritzenventil verbundenen Kapillaren und Schläuche werden mit der Waschflüssigkeit gefüllt und gespült.

#### Spülen der Spritze über Chromeleon

- 1. Öffnen Sie in Chromeleon das Dialogfenster Commands für den Autosampler.
- 2. Klicken Sie unter **Sampler** auf **PrimeSyringeNumber** und geben Sie ein, wie oft die Spritze zum Spülen gefüllt und entleert werden soll.
- 3. Geben Sie das Kommando **PrimeSyringe**, bis keine Blasen mehr sichtbar sind. Verwenden Sie gegebenenfalls frisch entgaste Waschflüssigkeit.
- 4. Wenn Sie die Luft auf diese Weise nicht aus der Spritze entfernen können, führen Sie die unter **Troubleshooting** genannten Schritte durch (siehe unten).

#### Spülen der Spritze vom Gerätedisplay aus

- 1. Wählen Sie die Funktionstaste **Prime**.
- 2. Entlüften Sie die Spritze, bis keine Blasen mehr sichtbar sind. Verwenden Sie gegebenenfalls frisch entgaste Waschflüssigkeit.
- 3. Wenn Sie die Luft auf diese Weise nicht aus der Spritze entfernen können, führen Sie die unter **Troubleshooting** genannten Schritte durch (siehe unten).

#### Troubleshooting

Wenn Sie die Luft nicht durch Spülen aus der Spritze entfernen können:

- 1. Bauen Sie die Spritze wie in Kapitel Wechseln der Spritze beschrieben aus  $(\rightarrow$  Seite 121).
- 2. Füllen Sie ein Becherglas mit der Waschflüssigkeit oder einer Mischung aus Isopropanol und Wasser (50:50).
- 3. Geben Sie die Spritze mit dem Spritzeneingang nach unten in das Becherglas.
- 4. Bewegen Sie die Spritze mit eingefahrenem Stößel mehrfach schnell hin und her. Damit werden Luftblasen entfernt, die sich unter Umständen am Spritzeneingang befinden.
- 5. Ziehen Sie die Spritze mehrfach hintereinander zügig auf und stoßen Sie den Inhalt schnell aus, bis die Spritze luftfrei ist.

- 6. Ziehen Sie die Spritze auf. Halten Sie die Spritze senkrecht und stoßen Sie etwas Flüssigkeit aus, damit sich am Spritzeneingang keine Luft mehr befindet.
- 7. Bauen Sie die Spritze wie in Wechseln der Spritze beschrieben wieder ein  $(\rightarrow$  Seite 121).
- 8. Führen Sie einen Waschzyklus durch. Führen Sie in Chromeleon den Befehl **Wash** aus oder wählen Sie am Gerätedisplay die Funktionstaste **Wash** oder wählen Sie **Wash** im Menü **Control** aus.

## 4.7 Einsetzen und Bestücken eines Probenhalters

Der Autosampler ist so gebaut, dass Sie die Probenhalter oder Mikrotiterplatten bequem einsetzen können. Eine Tabelle mit den verfügbaren Probenhaltern finden Sie auf Seite 23.



**Warnung:** Um Verletzungen zu vermeiden, greifen Sie während einer laufenden Analyse niemals in den Probenraum.

- 1. Schwenken Sie den Deckel des Drehtellers ganz nach links.
- 2. Holen Sie das Segment (rot, grün oder blau) nach vorn, in das Sie einen Probenhalter einsetzen möchten. Wählen Sie eine der folgenden Möglichkeiten:
  - Drehen Sie den Teller per Hand, bis das Segment vorn ist.
  - Wählen Sie die Funktionstaste Rotate, bis das Segment vorn ist.
  - Wählen Sie Rotate im Menü Control und wählen Sie dann R (G oder B).
  - Wählen Sie im Dialogfenster **Commands** in Chromeleon **InitiateChangeVial** und wählen Sie dann im Feld **Tray** den Eintrag **RedTray** (**GreenTray** oder **BlueTray**).
- Wenn Sie eine Mikrotiterplatte verwenden möchten Setzen Sie die Mikrotiterplatte auf dem Probenhalter an der Position A1 an und drücken Sie die Platte fest auf den Probenhalter, bis sie einrastet (→ Abb. 19).



Abb. 19: Aufsetzen einer Mikrotiterplatte auf einen Probenhalter

4. Setzen Sie den Probenhalter (oder den Probenhalter mit Mikrotiterplatte) in den Drehteller ein (→ Abb. 20). Der Probenhalter muss dabei auf den Positionierstiften einrasten.



Abb. 20: oben: Einsetzen und Bestücken eines Probenhalters unten: Drehteller (links Probenhalter mit Mikrotiterplatte, rechts mit Probenfläschchen)

An drei Positionen im Drehteller sind standardmäßig Halter für 5 Probenfläschchen à 10 mL installiert, z.B. zur Aufnahme der Reagensflüssigkeiten. Wenn Sie *ausnahms-weise ohne* diese Halter arbeiten möchten, achten Sie darauf, dass die Positionierklammer des Probenhalters rechts an der Farbmarkierung des jeweiligen Segments ausgerichtet ist.



Abb. 21: Ausrichtung des Probenhalters im Drehteller

- Bestücken Sie den Probenhalter gegebenenfalls mit den geeigneten Probenfläschchen (→ Seite 24).
- 6. Schwenken Sie den Deckel des Drehtellers wieder nach rechts, bis Kontakt zum Nadelarm besteht.

- 7. Wenn der Autosampler über Chromeleon gesteuert wird
  - a) Prüfen und ändern Sie in Chromeleon gegebenenfalls die Einstellungen für die Probengefäße (→ Seite 84). Die Probennadel kann beschädigt werden, wenn die Einstellungen in Chromeleon nicht mit den tatsächlichen installierten Probengefäßen übereinstimmen.
  - b) Passen Sie gegebenenfalls den Wert für die Nadelhöhe über den Parameter SampleHeight an (→ Seite 91). Beachten Sie die Informationen zu SampleHeightOffset auf Seite 93.

## 4.8 Bestücken der 5-Positionen Halter

An drei Positionen im Drehteller sind werkseitig Halter für jeweils 5 Probenfläschchen à 10 mL zur Aufnahme der Reagens- bzw. Transportflüssigkeiten installiert. Bestücken Sie den Halter mit den geeigneten 10 mL-Probenfläschchen ( $\rightarrow$  Seite 24).

Alternativ können Sie die Halter auch mit 2 mL-Probenfläschchen bestücken. In diesem Fall müssen Sie jedoch zuvor den Adapter (Best.-Nr. 6820.4092) installieren.



**Ing:** Um Verletzungen zu vermeiden, greifen Sie während einer laufenden Analyse niemals in den Probenraum.

- 1. Schwenken Sie den Deckel des Drehtellers ganz nach links.
- 2. Holen Sie das Segment (rot, grün oder blau), das Sie bestücken möchten, nach vorn. Wählen Sie eine der folgenden Möglichkeiten:
  - Drehen Sie den Teller per Hand, bis das Segment vorn ist.
  - Wählen Sie Funktionstaste Rotate, bis das Segment vorn ist.
  - Wählen Sie Rotate im Menü Control und wählen Sie dann R (G oder B).
  - Wählen Sie im Dialogfenster **Commands** in Chromeleon **InitiateChangeVial** und wählen Sie dann im Feld **Tray** den Eintrag **RedVials** (**GreenVials** oder **BlueVials**).
- 3. Bestücken Sie die Halter mit den Probenfläschchen (10 mL oder 2 mL) ( $\rightarrow$  Seite 24).
- 4. Schwenken Sie den Deckel des Drehtellers wieder ganz nach rechts, bis Kontakt zum Nadelarm besteht.
- 5. Wenn der Autosampler über Chromeleon gesteuert wird Prüfen und ändern Sie gegebenenfalls über den Parameter ReagentLiquidHeight den Wert für die Nadelhöhe (→ Seite 91). Beachten Sie insbesondere die Informationen zu SampleHeightOffset auf Seite 93. Bei Verwendung von 2-mL-Probenfläschchen muss der Wert für SampleHeightOffset\_T angepasst werden (→ Seite 92). Andernfalls wird die Probennadel beschädigt.

Verwenden Sie in allen fünf Positionen eines Halters Probenfläschchen desselben Typs.

## 4.9 Äquilibrieren des Systems

Äquilibrieren Sie das UltiMate 3000-System, *bevor* Sie den Autosampler zur Probenanalyse einsetzen:

- 1. Spülen Sie das gesamte System mit dem Anfangseluenten so lange, bis sich keine andere Flüssigkeit mehr im System befindet.
- 2. Heizen oder kühlen Sie alle temperaturgeregelten Module auf die für die Anwendung erforderliche Temperatur.
- 3. Stellen Sie am Detektor die Wellenlänge ein und schalten Sie die Lampen ein.
- 4. Beobachten Sie den Pumpendruck. Vergewissern Sie sich, dass der Druck für die jeweilige Anwendung korrekt ist und dass er stabil bleibt.
- 5. Beobachten Sie das Detektorsignal. Vergewissern Sie sich, dass Sie das für die Anwendung erwartete Basisliniensignal bekommen und dass das Signal stabil bleibt.

Führen Sie die Äquilibrierung über Chromeleon durch oder wählen Sie die für die Äquilibrierung erforderlichen Kommandos und Parameter über die Menüs der einzelnen Geräte aus.

#### Äquilibrieren des Systems über Chromeleon

- Wählen Sie die Befehle und Parameter im Dialogfenster Commands aus.
- Automatisieren Sie die Äquilibrierung, indem Sie ein Äquilibrierprogramm erstellen und ablaufen lassen (→ Seite 71).
- Verwenden Sie den SmartStartup-Assistenten (siehe unten), um das Äquilibrierprogramm zu erstellen und ablaufen zu lassen.

#### Erstellen eines Äquilibrierprogramms über den SmartStartup-Assistenten

- 1. Öffnen Sie den Assistenten über SmartStartup im Menü Batch.
- 2. Folgen Sie den Anweisungen auf den einzelnen Seiten des Assistenten. Klicken Sie auf die Schaltfläche **Hilfe**, wenn Sie weitere Informationen zu einer Seite benötigen.
- 3. Nach Beendigung des Assistenten
  - erstellt Chromeleon ein Äquilibierprogramm und die entsprechende Sequenz.
  - ♦ öffnet Chromeleon das Äquilibrierfenster f
    ür die auf der Zeitbasis installierten Ger
    äte (→ Abb. 22, Seite 61).
  - öffnet Chromeleon das Dialogfenster Start Batch on.

Klicken Sie auf Start, um mit der Äquilibrierung zu beginnen.

Das Äquilibrierfenster zeigt für jedes Gerät des Systems den Status der Äquilibrierung an.



Abb. 22: Äquilibrierfenster

#### Äquilibrieren des Systems über die Gerätemenüs

Wählen Sie die Befehle und Parameter in den Menüs der einzelnen Geräte aus. Informationen zu den Menüs des Autosamplers finden Sie im Kapitel 5.5.2 ( $\rightarrow$  Seite 75). Informationen zu den Menüs der anderen Systemmodule finden Sie in der *Bedienungsanleitung* des jeweiligen Gerätes.

# 5 Betrieb und Wartung

Der Autosampler kann über das Chromatographie-Management-System Chromeleon gesteuert werden. Informationen hierzu finden Sie im Kapitel 5.4 ( $\rightarrow$  Seite 67).

Zusätzlich stehen am Gerätedisplay Funktionstasten und Menüs zur Verfügung, über die bestimmte Aktionen direkt am Autosampler durchgeführt werden können, zum Beispiel während der Inbetriebnahme oder für Diagnose- oder Wartungsarbeiten. Sie können zum Beispiel:

- Bestimmter Funktionen ausführen (den Drehteller bewegen)
- Befehle ausführen (Waschen, Anfahren von Servicepositionen)
- Parameter einstellen (Leaksensormodus, Solltemperatur für die Probenthermostatisierung)
- Informationen zu Diagnosezwecken abfragen
- Gerätekonfiguration abfragen und ändern

Informationen hierzu finden Sie im Kapitel 5.5 ( $\rightarrow$  Seite 73).

## 5.1 Einschalten des Autosamplers

Schalten Sie den Autosampler zur Inbetriebnahme über den Netzschalter auf der Geräterückseite ein. Wenn der Autosampler über den Netzschalter eingeschaltet wird, passiert Folgendes:

- Auf dem Gerätedisplay erscheinen kurzzeitig allgemeine Informationen zum Autosampler: Gerätetyp, Seriennummer, Bootloader- und Firmware-Version.
- Der Autosampler führt einen Selbsttest durch, bei dem alle wesentlichen Baugruppen auf korrekte Funktion überprüft werden. Nach erfolgreichem Selbsttest erscheint die Statusanzeige auf dem Gerätedisplay (→ Seite 64).
- Tritt während des Selbsttests ein Fehler auf, ist der Autosampler nicht betriebsbereit. Die LED Status auf der Gerätevorderseite leuchtet rot und es erscheint eine Meldung auf dem Gerätedisplay. Wird der Autosampler unter Chromeleon betrieben, erscheint die Meldung auch im Chromeleon Audit Trail.
   Schalten Sie den Autosampler aus, korrigieren Sie den Fehler (→ Seite 103) und schalten Sie den Autosampler wieder ein.

Im Normalbetrieb brauchen Sie den Autosampler nicht über den Netzschalter auszuschalten. Verwenden Sie stattdessen die Standby-Taste auf der Gerätevorderseite (→ Seite 20). Drücken Sie die Taste ca. 1 Sekunde lang, damit der Autosampler den Modus ändert. Schalten Sie den Autosampler über den Netzschalter aus, wenn Sie dazu aufgefordert werden, zum Beispiel, für bestimmte Wartungsarbeiten.

## 5.2 Statusanzeige

Nach erfolgreichem Selbsttest erscheint die Statusanzeige auf dem Gerätedisplay:

Sample Position	Inject Volume
R-E3	0.100 μl
18 (10) °C	Loop: Inject

Abb. 23: Statusanzeige

Angezeigt werden

- Probenposition (Sample Position) Informationen zur Angabe der Probenposition finden auf Seite 65.
- Injektionsvolumen (Inject Volume)
- *Nur Autosampler mit Probenthermostatisierung* Aktuelle Temperatur und Solltemperatur (= Angabe in Klammern) für die Probenthermostatisierung
- Ventilposition (Loop: Inject oder Load)

Falls erforderlich, können Sie die Helligkeit der Hintergrundbeleuchtung und den Kontrast der Displayanzeige an Ihre Anforderungen anpassen ( $\rightarrow$  Seite 86).
# 5.3 Probenposition

Die einzelnen Segmente des Drehtellers sind durch die Farben Rot, Grün und Blau (R/G/B) gekennzeichnet. Diese Kennzeichnung ist bei der Angabe der Probenposition wichtig.



Abb. 24: Probenposition (hier: Rack-Preview aus Chromeleon)

Einzelheiten zur Angabe der Fläschchenpositionen in einem Probenhalter sowie zur Position der Kavitäten bei einer Mikrotiterplatte finden Sie unten. Informationen zur Positionsangabe der Fläschchen im 5-Positionen Halter finden Sie auf Seite 66.

**Hinweis:** Der Autosampler unterstützt die automatische Erkennung von Probenfläschchen und Mikrotiterplatten ( $\rightarrow$  Seite 95).

### Probenposition bei Probenfläschchen und Mikrotiterplatten

Bei der Probenposition steht der Buchstabe für das Segment, in dem sich das Probengefäß befindet an erster Stelle, also R, G oder B. Entsprechend einem Koordinatensystem folgen ein Buchstabe und eine Zahl. Dabei werden die einzelnen Reihen von innen nach außen mit einem Buchstaben bezeichnet. Die Zahl gibt die Position der Probe innerhalb einer Reihe an. Dabei erfolgt die Nummerierung von links nach rechts.



Rotes Segment (40 Vials)

Abb. 25: Beispiel Probennummerierung (hier: Probenhalter für 40 Probenfläschchen)

## Beispiel

Die Position R-C4 (in Chromeleon: RC4) gibt an, dass sich das Probenfläschchen an folgender Position befindet:

- rotes Segment (R)
- dritte Reihe (C)
- vierte Position von links (4)

## **Position im 5-Positionen Halter**

Bei der Angabe der Position für die 10-mL-Probenfläschchen steht der Buchstabe für das Segment, in dem sich das Probenfläschchen befindet, an erster Stelle (R, G oder B), gefolgt von einer Zahl von 1 bis 5.



Abb. 26: Positionen im 5-Positionen Halter

# Beispiel

Die Position B3 gibt an, dass sich das Probenfläschchen an folgender Position befindet:

- blaues Segment (B)
- dritte Position (3)

# 5.4 Steuerung über Chromeleon

Vergewissern Sie sich zunächst, dass folgende Voraussetzungen erfüllt sind:

- 1. Die Chromeleon-Software ist auf dem Rechner installiert und der Lizenzcode ist eingegeben.
- 2. Der Autosampler ist über eine USB-Verbindung mit dem Chromeleon-Rechner verbunden.
- 3. Der Autosampler ist in Chromeleon eingerichtet ( $\rightarrow$  Seite 37).

Damit der Autosampler über Chromeleon gesteuert werden kann, müssen Sie die Zeitbasis, auf welcher der Autosampler installiert ist, mit dem Chromeleon-Client verbinden  $(\rightarrow Kapitel 5.4.1)$ .

Die Steuerung kann auf zweierlei Art erfolgen:

- **Direkt** über die Parameter und Befehle im Dialogfenster **Commands** oder auf einem Steuerfenster (Control Panel) (→ Seite 68)
- Automatisch über ein Steuerprogramm (PGM) ( $\rightarrow$  Seite 71)

### 5.4.1 Verbinden mit Chromeleon

- 1. Starten Sie gegebenenfalls den Chromeleon Server Monitor und den Chromeleon Server ( $\rightarrow$  Seite 37).
- 2. Starten Sie den Chromeleon-Client über das Chromeleon-Symbol 🎏 auf dem Desktop.

Wenn das Chromeleon-Symbol nicht auf dem Desktop vorhanden ist, klicken Sie auf **Start** in der Taskleiste. Wählen Sie **Programme** (oder **Alle Programme**, abhängig vom Betriebssystem), wählen Sie dann **Chromeleon**, und klicken Sie danach auf **Chromeleon**.

3. Verbinden Sie den Chromeleon-Client mit der Zeitbasis, auf welcher der Autosampler installiert ist. Einzelheiten hierzu finden Sie für das Dialogfenster **Commands** auf der Seite 68 und für das Steuerfenster auf der Seite 69.

Wenn der Autosampler korrekt mit Chromeleon verbunden ist:

- leuchtet die LED Connected auf der Gerätevorderseite grün. •
- sind keine Eingaben über das Gerätedisplay möglich. Der Drehteller kann jedoch • weiterhin über die Funktionstaste Rotate bewegt werden.
- stehen weitere Funktionen zur Lebensdauerprognose von Verschleißteilen zur Verfügung •  $(\rightarrow$  Seite 88).
- stehen Diagnosetests zur Verfügung, mit denen der Autosampler und verschiedene seiner • Komponenten auf ihre Funktionsfähigkeit hin überprüft werden können ( $\rightarrow$  Seite 89)
- ist die Standby-Taste auf der Gerätevorderseite weiterhin aktiv. •

Trennen Sie den Autosampler immer über das Kommando Disconnect von Chromeleon, ehe Sie ihn am Netzschalter ausschalten.

## 5.4.2 Direkte Steuerung

Die Parameter und Befehle werden über das Dialogfenster Commands (F8-Box) eingegeben und ausgeführt. Direkte Befehle werden mit der Eingabe ausgeführt. Für den Routinebetrieb stehen die meisten Parameter und Befehle auch in einem Steuerfenster zur Verfügung.

#### Öffnen des Dialogfensters Commands für den Autosampler

- Öffnen Sie ein (beliebiges) Steuerfenster. Die Steuerfenster sind im Chromeleon-1. Browser in Verzeichnis Dionex Templates/Panels abgelegt und können mit einem Doppelklick geöffnet werden.
- 2. Verbinden Sie das Steuerfenster mit der Zeitbasis, auf welcher der Autosampler installiert ist. Wählen Sie im Menü Control den Befehl Connect to Timebase und legen Sie auf der Seite Timebase die Zeitbasis fest. Weitere Informationen zum Dialog Timebase erhalten Sie über die Schaltfläche Hilfe.



**i** Hinweis: Das Menü Control ist nur sichtbar, wenn ein Steuerfenster geöffnet ist

- 3. Drücken Sie F8 oder wählen Sie **Command** im Menü **Control**.
- Zeigen Sie die Parameter und Befehle für den Autosampler an, indem Sie auf das 4. Pluszeichen neben Sampler klicken.

Welche Parameter und Befehle angezeigt werden, hängt ab von

- ♦ der Chromeleon-Version
- den in der Autosamplerkonfiguration festgelegten Optionen ( $\rightarrow$  Seite 40).
- dem Anzeigefilter (Normal, Advanced oder Expert)

5. Ändern Sie den Anzeigefilter, falls erforderlich. Führen Sie in der Befehlsliste einen Rechtsklick aus und wählen Sie den Filter im Menü aus.



Abb. 27: Dialogfenster Commands

6. Vergewissern Sie sich, dass der Autosampler mit Chromeleon verbunden ist. Wenn er nicht verbunden ist, verbinden Sie den Autosampler über den Befehl **Connect**.

Eine Liste aller für den Autosampler verfügbaren Kommandos und Parameter finden Sie in der *Chromeleon-Hilfe*. Ergänzend zu den Autosamplerkommandos und Parametern können Sie über das Dialogfenster **Commands** auch auf alle Kommandos und Parameter der anderen Geräte zugreifen, die auf der ausgewählten Zeitbasis installiert sind.

### Öffnen des Steuerfensters für den Autosampler

1. Klicken Sie im Menü View auf Default Panel Tabset oder klicken Sie auf das entsprechende Symbol in der Werkzeugleiste 🔟 und verbinden Sie sich dann mit dem Chromeleon-Server.

Chromeleon erstellt zentrale Steuerfenster (Panel Tabsets;  $\rightarrow$  Abb. 28, Seite 70) für alle auf dem Server installierten Zeitbasen. Ein Panel Tabset enthält Steuerfenster für die einzelnen Geräte der Zeitbasis sowie ein oder mehrere Steuerfenster für systemweite Funktionen, z.B. für das Erstellen und Ausführung von Sequenzen. Weitergehende Informationen zu Panel Tabsets finden Sie in der *Chromeleon-Hilfe*.

- Klicken Sie auf dem Panel Tabset f
  ür Ihre Zeitbasis die Registerkarte Sampler an (→ Abb. 28).
- 3. Vergewissern Sie sich, dass der Autosampler mit Chromeleon verbunden (connected) ist (die LED der Schaltfläche ist grün). Wenn er nicht verbunden ist, verbinden Sie den Autosampler über die Schaltfläche **Connect**.



Abb. 28: Steuerfenster für den Autosampler innerhalb eines Panel Tabsets

Auf dem Control Panel stehen die Parameter und Befehle zur Verfügung, die für den Routinebetrieb des Autosamplers benötigt werden. Alle anderen Parameter und Befehle können Sie über das Dialogfenster **Commands** ausführen. Sie können das Dialogfenster direkt vom Panel Tabset aus über **Command** im Menü **Control** öffnen.

## 5.4.3 Automatische Steuerung

Beim automatischen Betrieb wird der Autosampler über ein von Ihnen erstelltes Programm (PGM) gesteuert. Das Programm können Sie automatisch mit einem Software-Assistenten erstellen oder manuell, indem Sie ein vorhandenes Programm editieren. Neben Programmen für die Probenanalyse können Sie auch Programme für andere Zwecke erstellen, zum Beispiel, benutzerdefinierte Programme für die Probenvorbereitung ( $\rightarrow$  Seite 96), Programme, um das HPLC-System automatisiert herunterzufahren ( $\rightarrow$  Seite 99), oder Programme, um sicherzustellen, dass das System nach einem Stromausfall wie gewünscht weiterarbeitet. Einzelheiten zu diesen Programmen finden Sie in der *Chromeleon-Hilfe*.

#### Erstellen eines Programms über den Programm-Assistenten

- 1. Rufen Sie den Programm-Assistenten auf. Wählen Sie dazu im Menü File den Befehl New und wählen Sie dann Program File aus der Liste.
- 2. Der Assistent führt Sie durch die Programmerstellung. Übernehmen oder ändern Sie gegebenenfalls auf jeder Seite des Assistenten die Einstellungen. Informationen zu den einzelnen Seiten erhalten Sie über die Schaltfläche Hilfe.
- 3. Nach Abschluss des Programm-Assistenten erstellt Chromeleon automatisch das Programm.
- 4. Starten Sie das Programm wie unten beschrieben ( $\rightarrow$  Seite 72).

#### **Manuelles Erstellen eines Programms**

1. Öffnen Sie ein vorhandenes Programm.

Öffnen Sie das Programm mit einem Doppelklick. - *oder* -

Wählen Sie im Menü File den Befehl **Open**. Wählen Sie in Dialogbox unter **Object of Type** den Eintrag **Program** und wählen Sie das Programm aus.



Abb. 29: Chromeleon Programm (hier: Programmansicht im Commands View)

2. Nehmen Sie in dem Programm die Änderungen vor.

Die Geräteansichten (Device Views) bieten die einfachste Möglichkeit, ein Programm zu ändern  $\rightarrow$  Abb. 29). Klicken Sie ein Gerät an und nehmen Sie die Änderungen auf den jeweiligen Geräteseiten vor. Die Eingaben werden direkt in Kommandos mit korrekter Syntax umgewandelt.

Wenn Sie einen Parameter in der Geräteansicht nicht einstellen können, können Sie in die Ansicht **Commands** wechseln und den Parameter dort editieren oder neu eingeben. Die Ansicht **Commands** zeigt das gesamte Programm mit den verschiedenen Befehlen in der zeitlichen Reihenfolge an. Weitere Informationen finden Sie in der *Chromeleon-Hilfe*.

3. Starten Sie das Programm wie unten beschrieben.

#### Starten eines Programms

#### Programm zur Probenanalyse

- 1. Erstellen Sie eine Probentabelle (Sequenz). Die Sequenz muss neben dem Programm auch eine Methode zur Auswertung der Probendaten enthalten (Peakidentifzierung, Flächen- und Stoffmengenbestimmung).
- 2. Weisen Sie das Programm und die Methode den einzelnen Proben in der Tabelle zu.
- 3. Nehmen Sie die Sequenz in den Batch auf und starten Sie den Batch.

Informationen zu den einzelnen Schritten finden Sie in der Chromeleon-Hilfe.

#### Andere Programme

Nehmen Sie die Programme in den Batch auf und starten Sie den Batch.

# 5.5 Funktionstasten und Menüs am Autosamplerdisplay

Über die Funktionstasten und Menüs des Gerätedisplays können Sie direkt am Autosampler verschiedene Einstellungen vornehmen und bestimmte Befehle ausführen.

Informationen zu den verschiedenen Funktionstasten finden Sie im Kapitel 5.5.1 (siehe unten) sowie auf der Seite 76. Informationen zu den einzelnen Menüs finden Sie im Kapitel 5.5.2 ( $\rightarrow$  Seite 75).

## 5.5.1 Einblenden der Funktionstasten

Vier weiße Punkte unterhalb des Displays markieren die Positionen der Funktionstasten **Menu**, **Rotate**, **Wash** und **Prime** ( $\rightarrow$  Tabelle auf Seite 74).

Berühren Sie den weißen Punkt ganz links auf der Glasplatte mit dem mitgelieferten Menüstift (Best.-Nr. 6300.0100), um die Funktionstasten einzublenden.



Abb. 30: Einblenden der Funktionstasten

Die Funktionstasten ersetzen die Informationen in der untersten Zeile der Statusanzeige. Wird keine Auswahl getroffen, erscheint nach circa 5 Sekunden wieder die ursprüngliche Zeile der Statusanzeige.



Abb. 31: Funktionstasten

Um	Wählen Sie
das Hauptmenü aufzurufen (→ Seite 77).	Menu
den Drehteller um ein Segment zu bewegen. Falls erforderlich, fährt die Probennadel zuvor aus dem Probenfläschchen (Mikrotiterplatte) heraus.	Rotate
einen Waschvorgang durchzuführen. Das Äußere der Probennadel wird im Waschport mit der Waschflüssigkeit gespült und danach über die Membranpumpe mit Luft getrocknet. Dabei wird das Volumen verwendet, dass Sie unter <b>Wash</b> im Menü <b>Control</b> festgelegt haben (→ Seite 78).	Wash
das Spritzenventil und die Spritze zu entlüften. Dabei wird die Spritze mehrfach aufgezogen und entleert. Weitere Informationen zum Spülen der Spritze finden Sie im Kapitel 4.6 ( $\rightarrow$ Seite 55).	Prime

Wenn der Autosampler in Chromeleon verbunden ist ("connected"), sind keine Eingaben über das Gerätedisplay möglich. Den Drehteller können Sie jedoch weiterhin über die Taste **Rotate** bewegen.

## 5.5.2 Autosampler-Menüs

Abb. 32 zeigt eine Übersicht über die verschiedenen Autosampler-Menüs. Informationen zum allgemeinen Aufbau der Menüs finden Sie auf Seite 76. Informationen zu den einzelnen Kommandos und Parametern, die in den verschiedenen Menüs zur Verfügung stehen, finden Sie in den Kapiteln 5.5.2.2 bis 5.5.2.6 ( $\rightarrow$  ab Seite 77).



Abb. 32: Menüstruktur (hier für Autosampler mit der Probenthermostatisierung)

# 5.5.2.1 Aufbau der Menüs

Die einzelnen Menüs sind wie folgt aufgebaut:



Abb. 33: Aufbau der Menüs (hier: Menü "Control")

Nr.	Beschreibung
1	Zeigt den Menünamen und die Anzahl der Menüpunkte an.
2	Die Menüpunkte werden als nummerierte Liste angezeigt. Der auswählbare Menüpunkt ist unterstrichen dargestellt.
3	Navigationsleiste

Je nach ausgewähltem Menüpunkt oder Parameter erscheinen unterschiedliche Tasten in der Navigationsleiste:

Um	Wählen Sie
zum vorherigen Eintrag in einer Liste zurückzugelangen. Sind mehr als 4 Punkte in der Liste vorhanden, können Sie nach Erreichen der 1. Zeile durch die Liste scrollen ( $\rightarrow$ Key autorepeat, Seite 80).	^
numerische Werte hochzuzählen.	^
zum nächsten Eintrag in einer Liste zu gelangen. Sind mehr als 4 Punkte in der Liste vorhanden, können Sie nach Erreichen der 4. Zeile durch die Liste scrollen ( $\rightarrow$ Key autorepeat, Seite 80).	V
zur nächsten Stelle einer Zahl weiterzugehen. Ein eventuell vorhandener Dezimal- punkt wird automatisch übersprungen.	>
die Auswahl zu bestätigen und gegebenenfalls das Eingabefeld zu aktivieren. Hat der Anwender nur Lesezugriff, ist die Funktionstaste <b>Select</b> nicht vorhanden	Select
eine Menüebene nach oben zu gelangen.	Back

Um	Wählen Sie
den Befehl auszuführen.	Execute
die Auswahl oder Eingabe zu bestätigen.	OK
die Aktion abzubrechen und den alten Wert wieder herzustellen.	Cancel
Hinweis: Abhängig vom ausgewählten Menüpunkt können spezifische Tasten die oben genannten Tasten in der Navigationsleiste ersetzen.	

Wird ein Fehler erkannt, blinken eine oder mehrere Meldungen auf dem Gerätedisplay. Dann erscheinen in der Navigationsleiste die Tasten **Prev**, **Next** und **Clear**.

Um	Wählen Sie
zur vorherigen Meldung zurück zu gelangen.	Prev
zur nächsten Meldung weiter zu gehen.	Next
die Meldung vom Gerätedisplay zu löschen.	Clear

### 5.5.2.2 Menü Main

Das Menü **Main** ist das Hauptmenü und damit die oberste Ebene in der Menüstruktur. Rufen Sie das Menü **Main** über die Funktionstaste **Menu** in der untersten Zeile der Statusanzeige ( $\rightarrow$  Seite 73) auf.

Über das Menü Main können Sie die folgenden Menüs aufrufen:

• Control

Im Menü **Control** können Sie Einstellungen für den Betrieb des Autosamplers vornehmen ( $\rightarrow$  Seite 78).

### • Maintenance

Im Menü **Maintenance** können Sie die für Wartungszwecke die erforderlichen Einstellungen vornehmen und Befehle ausführen ( $\rightarrow$  Seite 79).

### • Diagnostics

Im Menü **Diagnostics** erhalten Sie Informationen für Diagnosezwecke (Lesezugriff) und können einen Selbsttest für den Autosampler durchführen ( $\rightarrow$  Seite 79).

### • Configuration

Im Menü **Configuration** erhalten Sie Informationen zur Konfiguration des Autosamplers und können gegebenenfalls entsprechende Einstellungen vornehmen oder Befehle ausführen ( $\rightarrow$  Seite 80).

Der Autosamplertyp bestimmt, welche Kommandos und Parameter in den einzelnen Menüs zur Verfügung stehen.

# 5.5.2.3 Menü Control

Im Menü **Control** können Sie Einstellungen für den Betrieb des Autosamplers vornehmen. So können Sie das Volumen für den Waschvorgang festlegen, verschiedene Waschvorgänge starten und ein bestimmtes Segment im Drehteller nach vorn holen.

Um	Wählen Sie
das Volumen für den Waschvorgang festzulegen und einen Waschvorgang durchzuführen. Dabei wird das Äußere der Probennadel im Waschport mit Waschflüssigkeit gespült und nach dem Waschen über die Membranpumpe mit Luft getrocknet.	Wash
das Volumen für den Waschvorgang der Pufferschleife festzulegen und den Waschvorgang durchzuführen. Die Waschflüssigkeit wird über die Spritze aufgezogen, fließt durch die Pufferschleife und das Injektionsventil (Position Inject) und wird über die Injektor-Waste-Kapillare in den Abfall ausgestoßen.	Buffer loop wash
das Volumen für den Waschvorgang festzulegen und den Waschvorgang durchzuführen. Dabei wird das Innere der Probennadel mit Waschflüssigkeit gespült. Die Waschflüssigkeit wird über der Spritze aufgezogen, fließt durch die Pufferschleife, das Injektionsventil (Position Load) und die Probenschleife und wird durch die Probennadel in den Abfall ausgestoßen. Restflüssigkeit, die sich nach diesem Vorgang in der Probenschleife befindet, wird injiziert.	Inject loop wash
ein Segment auszuwählen (R = rot, G = grün, B = blau) und nach vorn zu holen. Der Probenhalter und der 5-Postionen Halter dieses Segments werden abwechselnd nach vorn geholt. Falls erforderlich, wird die Probennadel zuvor aus dem Probenfläschchen (Mikrotiterplatte) entfernt.	Rotate
die Temperatur für die Probenthermostatisierung festzulegen.	Nominal sample temp.
die Probenthermostatisierung ein- oder auszuschalten (On oder Off).	Sample cooling

## 5.5.2.4 Menü Maintenance

Über das Menü **Maintenance** können Sie für Wartungszwecke die erforderlichen Einstellungen vornehmen oder Befehle ausführen.

Um	Wählen Sie
die Spritze entsprechend der ausgewählten Option nach oben oder unten zu bewegen und das Volumen zu sehen, dass sich in der Spritze befindet.	Syringe
die Probennadel nach unten oder für Wartungszwecke in die Position <b>Maintain</b> zu bewegen.	Needle position
festzulegen, wohin die Waschflüssigkeit ausgestoßen wird.	Dispense destination
das Injektionsventil in die angegebene Position zu schalten (Load oder Inject).	Inject valve

## 5.5.2.5 Menü Diagnostics

Über das Menü **Diagnostics** erhalten Sie Informationen für Diagnosezwecke (Lesezugriff). Außerdem können Sie einen Selbsttest durchführen.

Um	Wählen Sie
einen Selbsttest durchzuführen. Wird ein Fehler gefunden, leuchtet die LED <b>Status</b> auf der Gerätevorderseite rot und eine Meldung erscheint auf dem Gerätedisplay.	Self test
die aktuelle Probentemperatur sehen.	Real sample temperature
die Temperatur des Kühlmittels zu sehen.	Coolant temperature
die Temperatur des Kühlkörpers zu sehen.	Heat sink temperature
die Kühlleistung zu sehen (in Prozent).	Cooling power
die Auslastung der Kühlpumpe zu sehen (in Prozent).	Coolant pump power
die Auslastung des Lüfters zu sehen (in Prozent).	Fan power
zu sehen, um welchen Autosampler es sich handelt.	Model
die installierte Firmware-Version zu sehen.	Firmware version
die Version des Bootloaders zu sehen.	Bootloader version
die Seriennummer des Autosamplers zu sehen.	Serial number

# 5.5.2.6 Menü Configuration

Über das Menü **Configuration** erhalten Sie Informationen zur Konfiguration des Autosamplers und können gegebenenfalls entsprechende Einstellungen vornehmen oder ändern.

Um	Wählen Sie
die Einstellungen für das Display und die Funktionstasten festzulegen: <b>Brightness</b> —legt die Helligkeit der Displayanzeige fest (in Prozent). <b>Contrast</b> —legt den Kontrast der Displayanzeige fest (in Prozent). <b>Key sound</b> —legt fest, ob bei Betätigung einer Funktionstaste ein akustisches Signal ertönt ( <b>On</b> = ja oder <b>Off</b> = nein). <b>Key autorepeat</b> — legt fest, ob bei längerer Tastenbetätigung der Tasten- druck automatisch wiederholt wird, z.B. zur schnellen Änderung eines Einstellwertes ( <b>On</b> = ja oder <b>Off</b> = nein).	Display & soft keys
<ul> <li>Festzulegen, ob und wie eine Leakerkennung erfolgen soll:</li> <li>Enabled—aktiviert die Leakerkennung. Wird eine Undichtigkeit erkannt, erscheint eine Meldung auf dem Gerätedisplay und es ertönt ein akustischer Alarm.</li> <li>Silent—aktiviert die Leakerkennung. Wird eine Undichtigkeit erkannt, erscheint nur eine Meldung auf dem Gerätedisplay. Es ertönt kein akustischer Alarm.</li> <li>Disabled—deaktiviert die Leakerkennung.</li> </ul>	Leak sensor mode
festzulegen, ob und wann bei der Injektion ein Waschvorgang durchgeführt werden soll:	Inject wash
<b>Before inject</b> —führt einen Waschvorgang aus, <i>bevor</i> die Probe aufgezogen wird (und damit <i>vor</i> der Injektion); die Probennadel enthält keine Probe. <b>After inject</b> —führt einen Waschvorgang aus, <i>nachdem</i> die Probe aufgezogen wurde (und bevor die Probennadel zum Injektionsport fährt); die Nadel wird von außen gewaschen.	
<b>Before &amp; after</b> —führt zwei Waschvorgänge durch: einen, <i>bevor</i> die Probe aufgezogen wird (die Probennadel enthält keine Probe) und einen, <i>nachdem</i> die Probe aufgezogen wurde (die Probennadel enthält Probe). <b>No wash</b> —führt keinen Waschvorgang aus.	
anzugeben, mit welchen Probengefäßen das jeweilige Segment bestückt ist.	Red (Green, Blue) tray
zu den Werkseinstellungen zurückzukehren. Es öffnet sich das Dialogfenster <b>Reset to factory defaults?</b> Bestätigen Sie diese Meldung mit <b>OK</b> , wenn Sie zu den Werkseinstellungen zurückkehren möchten. Brechen Sie die Aktion mit <b>Cancel</b> ab, wenn Sie Ihre Einstellungen beibehalten möchten.	Reset to factory defaults

# 5.6 Einstellungen für den Autosampler-Betrieb

In diesem Kapitel finden Sie Informationen zu Einstellungen und Funktionen, die Sie für den Betrieb des Autosamplers beachten sollten:

Informationen zum Betrieb	siehe Seite
Verbinden von Autosampler und Pumpe	Siehe unten
Synchronisieren von Injektionsbefehl und Kolbenhüben	82
Einschalten der Probentellerthermostatisierung	83
Ändern der Drehtellerkonfiguration	84
Festlegen der Injektionsparameter und Starten der Injektion	84
Erkennen von Undichtigkeiten	86
Anpassen von Helligkeit und Kontrast der Displayanzeige	86
SmartStartup und SmartShutdown	87
Empfehlungen für den Betrieb	87

Beachten Sie darüber hinaus auch die Informationen zu den speziellen Funktionen, die in Chromeleon zur Verfügung stehen ( $\rightarrow$  Seite 88).

# 5.6.1 Verbinden von Autosampler und Pumpe

Wenn das UltiMate 3000-System eine Pumpe aus der UltiMate 3000 Pumpenserie enthält, können Sie festlegen, mit welcher Pumpe der Autosampler verbunden werden soll. Nehmen Sie diese Einstellung grundsätzlich vor, da sie:

- angibt, welche Pumpe den Fluss liefert. Diese Information ist wichtig, wenn Sie im Bypass-Modus arbeiten (→ Seite 13).
- erlaubt, den Injektionsbefehl des Autosamplers mit den Kolbenhüben einer Niederdruckgradientenpumpe DPG-3600 bzw. LPG-3400 zu synchronisieren (→ Kapitel 5.6.2).

Im Programm **Server Configuration** legen Sie fest, mit welcher Pumpe der Autosampler standardmäßig verbunden ist:

- 1. Starten Sie das Programm Server Configuration ( $\rightarrow$  Seite 39).
- 2. Markieren Sie den Autosampler in der Zeitbasis mit einem Rechtsklick und wählen Sie im Menü den Punkt **Properties**.
- 3. Legen Sie auf der Registerkarte **Segments** / **Pump Link** unter **Pump Link** die Pumpe fest, mit der der Autosampler standardmäßig verbunden sein soll (→ Seite 42).

## 5.6.2 Synchronisieren von Injektionsbefehl und Kolbenhüben

Nur wenn das UltiMate 3000-System eine DGP-3600 oder LPG-3400 enthält

**i** Hinweis: Wenn es sich bei der Pumpe um eine LPG-3400XRS handelt *oder* wenn Sie eine LPG-3400XRS *und* eine weitere Pumpe mit dem Autosampler verbinden möchten, folgen Sie bitte der Beschreibung in der *Bedienungs-anleitung zur LPG-3400XRS*.

Nur wenn Sie den Autosampler im Programm **Server Configuration** mit einer DGP-3600 oder LPG-3400 verbunden haben ( $\rightarrow$  Kapitel 5.6.1), kann der Injektionszeitpunkt mit den Kolbenhüben der Pumpe synchronisiert werden. Damit wird sichergestellt, dass alle Injektionen zur selben Phase des Pumpenzyklus erfolgen. Auf diese Weise kann bei Gradientenanwendungen die Retentionszeitpräzision verbessert werden.

Im Dialogfenster **Commands**, im Programm (PGM) oder auf dem Steuerfenster für den Autosampler (unter **More Options**) können Sie festlegen, ob Sie mit oder ohne Synchronisierung arbeiten möchten und gegebenenfalls eine andere Pumpe festlegen:

- Wenn Sie ohne Synchronisierung arbeiten möchten, setzen Sie SyncWithPump auf Off.
- Unter **PumpDevice** wird der Name der Pumpe angezeigt, mit der die Synchronisierung erfolgt. Wenn die Synchronisierung mit einer anderen Pumpe erfolgen soll, geben Sie den Namen der Pumpe in das Eingabefeld ein. (Verwenden Sie den Namen, der im Programm **Server Configuration** für die Pumpe festgelegt ist.) Vergewissern Sie sich, dass **SyncWithPump** auf **On** gesetzt ist.

## 5.6.3 Einschalten der Probenthermostatisierung

#### Nur Autosampler mit Probenthermostatisierung

Die Probenthermostatisierung schalten Sie über Chromeleon oder am Gerätedisplay ein- und aus. Dort legen Sie auch die Temperatur fest.

#### Einschalten der Probenthermostatisierung über Chromeleon

- 1. Öffnen Sie das Dialogfenster Commands für den Autosampler.
- 2. Wählen Sie **Temperature** und **Nominal** und geben Sie die Temperatur ein. Durch die Eingabe einer Temperatur wird **TempCtrl** auf **On** gesetzt, falls noch nicht geschehen.



Abb. 34: Einschalten der Probenthermostatisierung

Wenn Sie für eine bestimmte Anwendung ohne Probenthermostatisierung arbeiten möchten, setzen Sie **TempCtrl** auf **Off**.

Wenn Sie die Probenthermostatisierung zu einem späteren Zeitpunkt wieder aktivieren möchten, setzen Sie **TempCtrl** auf **On** zurück. **TempCtrl** wird automatisch auf **On** gesetzt, wenn Sie unter **Nominal** eine andere Temperatur eingeben.

#### Einschalten der Probenthermostatisierung am Gerätedisplay

- 1. Rufen Sie am Gerätedisplay das Menü Control auf.
- 2. Wählen Sie Nominal sample temperature und geben Sie die Temperatur ein.
- 3. Wählen Sie Sample Cooling und On, um die Probenthermostatisierung einzuschalten.

Setzen Sie **Sample Cooling** auf **Off**, wenn Sie zu einem späteren Zeitpunkt ohne Probenthermostatisierung arbeiten möchten.

# 5.6.4 Ändern der Drehtellerkonfiguration

Bei der Installation des Autosamplers in Chromeleon wird im Programm **Server Configuration** auf der Seite **Segments / Pump Link** festgelegt, welche Probengefäße standardmäßig im Drehteller installiert sind. Wenn Sie zu einem späteren Zeitpunkt andere Probengefäße im Autosampler installieren, müssen Sie die Einstellung in Chromeleon entsprechend anpassen:

- Wenn Sie die Standardeinstellung ändern möchten Nehmen Sie die Änderung im Programm Server Configuration auf der Seite Segments/ Pump Link vor (→ Seite 41).
- *Wenn Sie die Standardeinstellung nicht ändern möchten* Sie können diese Anpassung direkt im Chromeleon Client vornehmen. Wählen Sie eine der folgenden Alternativen:
  - Öffnen Sie das Dialogfenster **Commands** und legen Sie unter **RedTray**, **GreenTray** und **BlueTray** fest, mit welchen Probengefäßen das jeweilige Segment bestückt ist.
  - Öffnen Sie das Steuerfenster für den Autosampler und legen Sie die Probengefäße über **Tray Type** unter **Red Section**, **Green Section** und **Blue Section** fest.

Stimmen die Konfiguration des Autosamplers und die Einstellungen in Chromeleon nicht überein, kann es zu einer Beschädigung der Probennadel kommen.

### 5.6.5 Festlegen der Injektionsparameter und Starten der Injektion

In Chromeleon können Sie den Injektionsbefehl manuell geben oder in ein Programm aufnehmen. Wenn die Injektion erfolgt ist, sendet der Autosampler sendet ein Signal an Chromeleon.

Die allgemeinen Injektionsparameter sind in Chromeleon voreingestellt und für Wasser als Lösungsmittel optimiert, z.B. **DrawSpeed** (= Geschwindigkeit, mit der die Probe aufgezogen wird), **DispSpeed** (= Geschwindigkeit, mit der die Spritze über die Pufferschleife ausstößt), und **DrawDelay** (= Zeit, die die Probennadel nach dem Aufziehen der Probe noch im Probenfläschchen bleibt). Prüfen und ändern Sie diese Parameter entsprechend Ihrer Anwendung ab.

## Hinweis: Wenn die Probe, der Eluent und/oder die Reinigungslösung einen hohen Gasgehalt aufweisen, wählen Sie eine niedrigere DrawSpeed und einen längeren DrawDelay.

#### Injektionsbefehl manuell geben

**i** Hinweis: Vergewissern Sie sich, dass als Injektionsmodus Normal ausgewählt ist. Prüfen und ändern können Sie dies im Dialogfenster Commands unter Inject Mode oder auf dem Steuerfenster für den Autosampler (über More Options unter General/Injection).

Um die Injektionsparameter einzugeben und die Injektion zu starten, wählen Sie eine der folgenden Alternativen,

- Dialogfenster Commands
   Wählen Sie Inject. Legen Sie unter Position die Probenposition fest. Legen Sie unter Volume das Injektionsvolumen fest. Starten Sie die Injektion mit Execute.
- Steuerfenster für den Autosampler Legen Sie im Abschnitt **Inject** die Probenposition und das Injektionsvolumen fest. Starten Sie die Injektion mit **Inject**.
- Menü Control oder Symbolleiste Online

Das Menü **Control** und die Symbolleiste **Online** sind nur verfügbar, wenn ein Steuerfenster geöffnet und mit der Zeitbasis verbunden ist, auf welcher der Autosampler installiert ist.

Wählen Sie **Inject** im Menü **Control** oder klicken Sie auf das Symbol **Inject** auf der Symbolleiste **Online**. Legen Sie im Dialogfenster die Probenposition und das Injektionsvolumen fest. Starten Sie die Injektion mit **OK**.

#### Injektionsbefehl über ein Programm geben

Damit die Injektion ausgeführt werden kann, muss das Programm folgende Zeile enthalten 0.000 Inject.

Diese Zeile besagt, dass der Injektionsbefehl zur Retentionszeit 0.000 Minuten ausgeführt wird. Wenn Sie das Programm über den Programm-Assistenten erstellen ( $\rightarrow$  Seite 71), wird dieser Injektionsbefehl automatisch in das Programm aufgenommen.

Legen Sie die Probenposition und das Injektionsvolumen in der Sequenz fest und starten Sie das Programm ( $\rightarrow$  Seite 72).

# 5.6.6 Erkennen von Undichtigkeiten (Leakerkennung)

Die Leakerkennung aktivieren und deaktivieren Sie in Chromeleon oder am Gerätedisplay. Wird eine Undichtigkeit erkannt, führt dies *nicht* zum Abbruch der Analyse.

#### Einschalten der Leakerkennung über Chromeleon

- 1. Öffnen Sie das Dialogfenster Commands für den Autosampler.
- 2. Wählen Sie LeakSensorMode und dann Enabled oder Silent.

Enabled-aktiviert die Leakerkennung. Wird eine Undichtigkeit erkannt

- leuchtet die LED Status auf der Gerätvorderseite rot.
- erscheint eine Meldung in Chromeleon und auf dem Gerätedisplay.
- ertönt ein akustischer Alarm.

Silent—aktiviert die Leakerkennung. Wird eine Undichtigkeit erkannt

- leuchtet die LED **Status** auf der Gerätvorderseite rot.
- erscheint eine Meldung in Chromeleon und auf dem Gerätedisplay.
- ertönt jedoch *kein* akustischer Alarm.

Über Disabled wird die Leakerkennung ausgeschaltet.

#### Einschalten der Leakerkennung am Gerätedisplay

- 1. Wählen Sie die Funktionstaste Menu.
- 2. Rufen Sie das Menü Configuration auf und wählen Sie Leak sensor mode.
- 3. Wählen Sie Enabled, Silent oder Disabled (siehe oben).

### 5.6.7 Anpassen von Helligkeit und Kontrast der Displayanzeige

Die Helligkeit der Hintergrundbeleuchtung und den Kontrast der Displayanzeige können Sie in Chromeleon oder am Gerätedisplay an Ihre Anforderungen anpassen:

- Öffnen Sie in Chromeleon das Dialogfenster Commands für den Autosampler
   (→ Seite 68). Ändern Sie unter Brightness den Wert für die Helligkeit der Hintergrundbeleuchtung; ändern Sie unter Contrast den Wert für den Kontrast der Displayanzeige.
- Rufen Sie am Gerätedisplay das Menü **Configuration** und dann **Display & soft keys** auf (→ Seite 80). Ändern Sie unter **Brightness** den Wert für die Helligkeit der Hintergrundbeleuchtung; ändern Sie unter **Contrast** den Wert für den Kontrast der Displayanzeige.

# 5.6.8 SmartStartup und SmartShutdown

Der SmartStartup-Assistent hilft Ihnen, wiederkehrende Tätigkeiten zu automatisieren ( $\rightarrow$  Seite 60). SmartStartup übernimmt das automatisierte und kontrollierte Einschalten der verschiedenen Module Ihres UltiMate 3000-Systems. So werden beispielsweise beim Autosampler mit Probenthermostatisierung der Drehteller und die darin befindlichen Komponenten auf die eingestellte Temperatur gebracht. Zusätzlich kann SmartStartup die Pumpe automatisch entlüften, die Trennsäule spülen und das HPLC-System Äquilibrieren. Wichtige Modulparameter werden automatisch kontrolliert. Wenn die Grenzwerte von den Modulen eingehalten werden, kann die zuvor aufgesetzte Probensequenz automatisch gestartet werden. SmartStartup kann zu einem beliebigen Zeitpunkt aktiviert werden.

Mit dem **SmartShutdown**-Assistenten können Sie ein Programm erstellen, mit dem das HPLC-System zur kurzfristigen Außerbetriebnahme in den Bereitschafts-Modus (Standby) versetzt oder für eine längere Betriebsunterbrechung automatisiert heruntergefahren (Shutdown) werden kann ( $\rightarrow$  Seite 99).

# 5.6.9 Empfehlungen für den Betrieb

Führen Sie folgende Aktionen in regelmäßigen Abständen durch, besonders nach längeren Stillstandszeiten:

- Spülen Sie die Spritze ( $\rightarrow$  Seite 55).
- Führen Sie einen Waschzyklus für die Pufferschleife durch. Verwenden Sie in Chromeleon den Befehl WashBufferLoop und am Gerätedisplay im Menü Control der Befehl Buffer loop wash. Verwenden Sie zum Waschen der Pufferschleife mindestens das doppelte Volumen der Pufferschleife. Eventuell müssen Sie die Spritze öfters aufziehen.

# 5.7 Spezielle Funktionen in Chromeleon

Dieses Kapitel gibt einen kurzen Überblick über einige spezielle Funktionen, die für den Autosampler in Chromeleon zur Verfügung stehen.

Erfahren Sie mehr über	Auf Seite
Überwachung von Verschleißteilen	siehe unten
Diagnosetests	90
Definieren der Nadelhöhe (Sample Height, Reagent Liquid Height)	91
Bewegen des Drehtellers vor dem Aufziehen der Probe (Tray Shake)	94
Belüften des Probenfläschchens (Puncturer Offset)	94
Operational Qualification und Performance Qualification	94
Automatischer Tray Test	94
Automatische Probenfläschenerkennung	95
Benutzerdefinierte Programme	96

Diese Funktionen können Sie (soweit nicht anders angegeben) über das Dialogfenster **Commands** aufrufen. Zusätzlich stehen einige dieser Funktionen auch auf dem Steuerfenster für den Autosampler zur Verfügung. Weitergehende Informationen zu den genannten Funktionen finden Sie in der *Chromeleon-Hilfe*.

# 5.7.1 Aktive Überwachung von Verschleißteilen (Predictive Performance)

Predictive Performance (= aktive Überwachung der wichtigsten Verschleißteile) unterstützt Funktionen zur Lebensdauerprognose von Verschleißteilen sowie zur Kontrolle und Dokumentation von Service- und (Re-)Qualifizierungsmaßnahmen.

#### **Dialogfenster Commands**

Öffnen Sie das Dialogfenster **Commands** für den Autosampler und legen Sie die Grenzwerte fest. Eine vollständige Liste der für den Autosampler verfügbaren Befehle und Zähler finden Sie in der *Chromeleon-Hilfe*. Um die Informationen für die Predictive Performance aktuell zu halten, führen Sie folgende Befehle aus ( $\rightarrow$  Tabelle).

Geben Sie nach	folgenden Befehl
einem Tausch des Nadelsitzes	NeedleSeatChanged
einem Tausch der Probennadel	NeedleChanged
einem Tausch der Rotordichtung	RotorSealChanged

Geben Sie nach	folgenden Befehl
einem Tausch der Spritze	SyringeChanged
einem Service (z.B. jährliche Wartung)	ServiceDone
einer Qualifizierung	QualificationDone

Über diese Befehle werden die Zähler zurückgesetzt und das Datum eingetragen, an dem die Maßnahme erfolgt ist.

#### Steuerfenster

Auf dem Steuerfenster für den Autosampler stehen Befehle und Zähler für die Predictive Performance über die Schaltflächen **Wellness**, **Qualification** und **Service** zur Verfügung. Hier können Sie die Grenzwerte eingeben und die Zähler zurücksetzen. Darüber hinaus zeigen Statusbalken die Qualifizierungs- und Serviceintervalle optisch an (ab Chromeleon 6.80).

Die Farbkodierung der Balken gibt Auskunft über den jeweiligen Status:

Farbe	Beschreibung	
Grün	OK.	
Gelb	Der Grenzwert ist fast erreicht oder das entsprechende Teil sollte demnächst gewartet oder getauscht werden.	
Orange	(Nur bei Anzeigen für die Eigenschaft "Qualification"). Der Grenzwert ist erreicht. Es gibt jedoch noch eine Toleranzfrist (Grace Period), in der das Gerät weiter verwendet werden darf.	
Rot	Der Grenzwert ist erreicht (bei Qualification: das Ende der Grace Period) und ein Austausch, Service oder Qualifizierung des Gerätes ist überfällig. Das Gerät kann nicht mehr betrieben werden. Darüber hinaus ist es auch nicht möglich, einen Batch zu starten.	

Wird ein Grenzwert erreicht, erscheint außerdem eine Meldung im Chromeleon Audit Trail.

# 5.7.2 Diagnosetests

In Chromeleon stehen Diagnosetests zur Verfügung, mit denen Sie den Autosampler und verschiedene seiner Komponenten auf ihre Funktionsfähigkeit hin prüfen können. Das Diagnosetool-Kit für die UltiMate 3000-Autosampler (Best.-Nr. 6822.0030) enthält alle Materialien, die für die Durchführung der Tests benötigt werden.

- 1. Wählen Sie **Diagnostics** im Menü **Control**. (Das Menü **Control** ist nur sichtbar, wenn ein Steuerfenster geöffnet ist.)
- 2. Im Dialogfenster **Diagnostics** werden die Diagnosetests für alle Geräte angezeigt, die auf der aktuellen Zeitbasis installiert sind. Wählen Sie einen Test für den Autosampler aus. Hinweise zur Durchführung der Tests finden Sie in der *Chromeleon-Hilfe*.

Prüfen Sie	Mit folgendem Test
die Schraubverbindungen sowie den Autosampler insgesamt auf Dichtigkeit. Der Test kann nur in Verbindung mit einer Pumpe des UltiMate 3000-Systems durchgeführt werden.	Leak Test
alle Antriebe und Sensoren des Autosamplers auf korrekte Funktionsweise.	Drives and Sensors Test

Wurde ein Test nicht bestanden, finden Sie im Kapitel Diagnose-Meldungen Informationen zu möglichen Ursachen sowie Vorschläge für Abhilfemaßnahmen ( $\rightarrow$  Seite 110).

## 5.7.3 Definieren der Nadelhöhe (Sample Height, Reagent Liquid Height)

Der Parameter **SampleHeight** beschreibt den Abstand (in mm) zwischen dem inneren Boden des Probengefäßes (Probenfläschchen, Mikrotiterplatte, Deep-Well-Plate) und der Nadelspitze. Dieser Wert definiert, wie tief die Probennadel beim Ansaugen der Probe in das Probengefäß eintaucht.

Der Parameter **ReagentLiquidHeight** definiert diesen Wert, für die 10-mL-Probenfläschchen im 5-Positionen Halter.



Nr.	Beschreibung
1	Probennadel
2	Probengefäß
3	Höhe der Probennadel über dem inneren Boden des Probengefäßes (SampleHeight/ReagentLiquidHeight)
4	Ausgleichswert für Bodenhöhe (SampleHeightOffset_XX)
5	Probenhalter
6	Drehteller

Abb. 35: Definieren der Nadelhöhe a) für Standard-Probenfläschchen b) für Standard-Mikrotiterplatte c) für Deep-Well-Plate

Bei **SampleHeight** (**ReagentLiquidHeight**) = **0** erreicht die Probennadel den Boden des Probengefäßes. Verschiedene Probengefäßtypen können jedoch eine unterschiedliche absolute Bodenhöhe haben. Daher könnte die Probennadel unbeabsichtigt auf den Boden des Probengefäßes auftreffen, wenn Sie zwischen verschiedenen Probengefäßtypen wechseln. Dabei kann die Probennadel beschädigt werden. Die Nadelspitze ist sehr empfindlich und kann sich leicht verformen.

Damit Sie für alle unterstützten Probengefäßtypen die gleiche **SampleHeight**-Einstellung verwenden können, steht für jeden Probengefäßtyp ein Wert zum Ausgleich der Bodenhöhe (= **SampleHeightOffset\_XX**) zur Verfügung (→ Tabellen auf Seite 92). Wenn die Probennadel in einen bestimmten Typ Probengefäß eintaucht, berücksichtigt Chromeleon diesen Wert entsprechend der Drehtellerkonfiguration. Mit den voreingestellten Werten berührt die Probennadel den Boden des Probengefäßes auch dann nicht, wenn SampleHeight (ReagentLiquidHeight) auf 0 mm gesetzt ist.

Bei Auslieferung des Autosamplers sind die in der Tabelle genannten Werte voreingestellt:

Für Probengefäß	SampleHeightOffset_XX [mm]
1,2-mL-Probenfläschchen	SampleHeightOffset_72 = 2.0
2-mL-Probenfläschchen	SampleHeightOffset_ $40 = 2.0$
4-mL-Probenfläschchen	SampleHeightOffset_22 = 2.0
10-mL-Probenfläschchen (im Probenhalter)	SampleHeightOffset_10 = 2.0
10-mL-Probenfläschchen (im 5-Positionen Halter)	SampleHeightOffset_T = $2.0$
Mikrotiterplatte (12 - 24 mm hoch) mit 96 Kavitäten	SampleHeightOffset_96 = 4.0
Mikrotiterplatte (12 - 24 mm hoch) mit 384 Kavitäten	SampleHeightOffset_384 = 3.0
Deep-Well-Plate (34 - 46 mm hoch) mit 24 Kavitäten	SampleHeightOffset_24deep = 2.0
Deep-Well-Plate (34 - 46 mm hoch) mit 96 Kavitäten	SampleHeightOffset_96deep = 2.0
Deep-Well-Plate (34 - 46 mm hoch) mit 384 Kavitäten	SampleHeightOffset_384deep = 2.0

Bei Verwendung der folgenden Probengefäße müssen Sie die Werte anpassen (→ Tabelle):

Für Probengefäß	SampleHeightOffset_XX [mm]
0,3- mL-Mikrodialyse-Probenfläschchen	SampleHeightOffset_72 = 13.0
0,3- mL zylindrische Probenfläschchen	SampleHeightOffset_72 = 2.0
0,5-mL-Probenfläschen (Eppendorf)	SampleHeightOffset_40 = 14.0
1,5-mL-Probenfläschen (Eppendorf)	SampleHeightOffset_40 = 2.0
2-mL-Probenfläschchen (bei Verwendung mit Adapter 6820. 4092 im 5-Positionen Halter)	SampleHeightOffset_T = $9.0$
Deep-Well-Plate (30 - 36 mm hoch) mit 96 Kavitäten	SampleHeightOffset_96deep = 7.0
Deep-Well-Plate (30 - 36 mm hoch) mit 384 Kavitäten	SampleHeightOffset_384deep = 7.0
Deep-Well-Plate (20 - 32 mm hoch) mit 96 Kavitäten	SampleHeightOffset_96deep = 12.0
Deep-Well-Plate (20 - 32 mm hoch) mit 384 Kavitäten	SampleHeightOffset_384deep = 12.0
Low-Well-PCR-Plate mit 384 Kavitäten (8 - 12 mm hoch)	SampleHeightOffset_384 = 8.0

⚠ Vorsicht:	Um eine Beschädigung der Probennadel zu vermeiden, dürfen die genannten Werte für <b>SampleHeightOffset_XX</b> keinesfalls unter- schritten werden. Dies gilt besonders, wenn Sie für <b>SampleHeight</b> ( <b>ReagentLiquidHeight</b> ) einen sehr kleinen Wert oder <b>0</b> gewählt haben.
<b>i</b> Hinweis:	Wenn Sie aus einem Probengefäß nur <i>ein</i> Mal injizieren möchten, wählen Sie den Wert für <b>SampleHeight</b> so, dass die Probennadel nur so tief wie nötig in das Probengefäß eintaucht. Damit wird eine Ver- schleppung zwischen den Proben weiter reduziert.

# 5.7.4 Bewegen des Drehtellers vor dem Aufziehen der Probe (Tray Shake)

Zur Rehomogenisierung der Probe, speziell nach längeren Standzeiten, unterstützt Chromeleon den "Tray Shake". Durch schnelle Drehtellerbewegungen wird die Probe vor dem Aufziehen gleichmäßig durchmischt. Tray Shake verhindert Peakflächenänderungen durch eine lokale Änderung der Konzentration im Probengefäß (z.B. durch Kondensation im Headspace).

Die Funktion ist standardmäßig deaktiviert (**AutoTrayShakeTimes** = 0). Geben Sie einen Wert ein um festzulegen, wie viele Schüttelbewegungen vor dem Aufziehen der Probe ausgeführt werden sollen.

## 5.7.5 Belüften des Probenfläschchens

Zur Verbesserung der Reproduzierbarkeit geben Sie einen Wert für den Parameter **PunctureOffset** ein.

Der Drehteller wird um den Wert verdreht, während sich die Probennadel im Septum befindet. Das Einstichloch wird entsprechend größer, so dass ein Druckausgleich stattfindet. Bei Probenfläschchen mit ungeschlitzten Septen und bei großen Injektionsvolumina ist dies besonders wichtig.

# 5.7.6 Operational Qualification und Performance Qualification

Um die Leistungsfähigkeit des HPLC-Systems zu kontrollieren und dokumentieren, führen Sie die Operational und Performance Qualification durch. Alle erforderlichen Materialien und eine detaillierte Anleitung zur Durchführung sind auf Anfrage erhältlich.

# 5.7.7 Automatischer Tray-Test

Der automatische Tray-Test prüft, ob die tatsächliche Bestückung des Probenhalters, aus dem die nächste Injektion erfolgen soll, der festgelegten Konfiguration entspricht. Ist dies nicht der Fall, kann die Probennadel beschädigt werden. Der Tray-Test wird vor jeder ersten Injektion aus einem Probenhalter nach einem manuellen Eingriff einmal pro Probenhalter ausgeführt.

Der Tray-Test ist standardmäßig aktiviert. Wenn Sie diese Funktion nicht nutzen möchten, setzen Sie das Property **EnableTrayTest** in Chromeleon auf **Disabled**. Ist der Niederhalter-Adapter (Best.-Nr. 6820.2402) installiert, *müssen* Sie den automatischen Tray-Test ausschalten.

# 5.7.8 Automatische Probengefäßerkennung

Ein Sensor im Niederhalter ( $\rightarrow$  Abb. 3, Seite 19) erkennt automatisch Probenfläschchen, Mikrotiterplatten und Deep-Well-Plates.

Der Sensor prüft, ob der Niederhalter auf einem Probengefäß aufsetzt. Befindet sich an einer angegebenen Position kein Probengefäß, erscheint eine Meldung am Gerätedisplay und im Chromeleon Audit Trail.

## 5.7.9 Benutzerdefinierte Programme

#### 5.7.9.1 Benutzerdefinierte Programme erstellen

Mit dem Programm-Assistenten von Chromeleon können Sie auch benutzerdefinierte Programme erstellen, zum Beispiel, um die einzelnen Schritte zur Probenvorbereitung festzulegen.

- 1. Rufen Sie den Programm-Assistenten auf ( $\rightarrow$  Seite 71).
- 2. Klicken Sie auf der Seite Sampler Options unter Inject Mode auf UserProg.

Program Wizard: Sampler Options			
	Inject <u>M</u> ode:	Nomal	•
	Connected Pump Device	Nomal	
	Bypass	L <u>S</u> ynchronize injec	tion with pump
	Flush Out Factor		[1,010,0]
	Maximum <u>I</u> nject Volume		[0,001100,000 µl]
	Flow		
	Bypass Time		[min]

Abb. 36: Benutzerdefiniertes Programm

3. Legen Sie auf der nächsten Seite des Assistenten die Schritte zur Probenvorbereitung fest.

Der Befehl **Inject** im Programm zur Probenanalyse startet das benutzerdefinierte Programm. Damit wird der chromatographische Lauf angehalten, bis das benutzerdefinierte Programm der Befehl **UDPInjectMarker** sendet.

Weitere Informationen und Beispiele für benutzerdefinierte Programme finden Sie in der *Chromeleon-Hilfe*.

**i** Hinweis: Die Befehle für die Probenvorbereitung (Udp-Kommandos) stehen auch im Dialogfenster **Commands** zur Verfügung.

# 5.7.9.2 Eingabe relativer Probenpositionen

In benutzerdefinierten Programmen können Sie Probenpositionen auch relativ zu einer anderen Position angeben. Beispielweise können Sie eine relative Position für das Probenfläschchen angeben, das zur Probenvorbereitung verwendet werden soll (PrepVial).

Die relative Position kann sich in einer anderen Reihe auf dem Probenhalter, in einem anderen Drehtellersegment, oder eine bestimmte Anzahl von Positionen von einem anderen Probenfläschchen entfernt befinden.

Informationen, wie Sie relative Probenpositionen in einem Programm eingeben müssen, finden Sie in der *Chromeleon-Hilfe*.

# 5.8 Außerbetriebnehmen des Autosamplers

Beachten Sie die folgenden Hinweise für die Außerbetriebnahme und den Transport des Autosamplers:

- Falls erforderlich, spülen Sie den Autosampler von Lösungsmitteln frei. Wenn ein Puffer als Teil der mobilen Phase eingesetzt wird, spülen Sie das System vor der Außerbetriebnahme einige Male mit einem Gemisch aus Methanol und Wasser (50:50). Dadurch vermeiden Sie eine Aufkonzentrierung von Salzen im System.
- Bei Betriebsunterbrechungen von 1 Woche oder länger, füllen Sie den Autosampler mit Methanol oder einem ähnlichen Alkohol wie Isopropanol oder Ethanol. Sind die Lösungsmittel im Autosampler nicht wasserlöslich, müssen Sie diese schrittweise ersetzen.
- Um den Autosampler zu transportieren oder zu versenden, müssen Sie
  - die Drainage entleeren. Schalten die Membranpumpe mit dem Befehl **PurgePumpOn** in Chromeleon für eine kurze Zeit ein.
  - die beweglichen Teile des Drehtellers (Probenfläschchen, Probenhalter, Mikrotiterplatten) und die Flasche für die Waschflüssigkeit entnehmen.
  - ♦ den Nadelarm und den Drehtellerdeckel beschrieben über das Schaumteil sichern (→ Verpackungsvorschrift).
  - ▲ Vorsicht: Dies ist ein Muss! Um eine Beschädigung der Probennadel und des Waschports zu vermeiden, fahren Sie die Probennadel und den Nadelarm in die korrekte Position (Nadel rechts und Nadel nicht im Nadelsitz oder Waschport). Drücken Sie die Taste Standby auf der Gerätevorderseite oder geben Sie den Befehl Standby in Chromeleon. Schalten Sie den Autosampler aus und sichern Sie den Nadelarm über das Verpackungsschaumteil.
- Versenden Sie das Gerät immer in der Originalverpackung. Beachten Sie die Verpackungsvorschrift.

Ist die Originalverpackung nicht mehr verfügbar, können Sie geeignete Geräteverpackungen über die Thermo Fisher Scientific-Vertriebsorganisation für Dionex HPLC-Produkte bestellen. Die Verpackungsvorschrift ist im Ordner "Installation and Qualification Documents for Chromatography Instruments" enthalten. Sie ist ebenfalls auf Anfrage erhältlich.

Erfolgt der Versand des Gerätes nicht in der Originalverpackung, entfällt die Gerätegarantie.

Wenn der Autosampler über Chromeleon gesteuert wird, können Sie den Autosampler und das HPLC-System über Chromeleon in den Standby-Modus setzen oder automatisiert herunterfahren ( $\rightarrow$  Seite 99).

#### Standby-Programm

Ein Standby-Programm versetzt das HPLC-System in den Bereitschaftszustand. Die wichtigsten Programmschritte sind:

- Am Programmende verringert das Programm automatisch den Fluss.
- Es fährt die Temperatur aller temperaturgesteuerten Systemmodule herunter.

Aus dem Bereitschaftszustand heraus können Sie die Applikationen schnell wieder reaktivieren.

#### Shutdown-Programm

Ein Shutdown-Programm fährt das HPLC-System automatisch herunter. Die wichtigsten Programmschritte sind:

- Das Programm verringert automatisch den Fluss am Programmende.
- Es schaltet bestimmte Komponenten und Funktionen des Systems aus (z.B. Detektorlampen, Temperaturregelung).

#### Erstellen eines Standby- oder Shutdown-Programms

Wählen Sie eine der folgenden Alternativen:

- Wählen Sie die Befehle und Parameter im Dialogfenster Commands aus.
- Automatisieren Sie die Außerbetriebnahme, indem Sie ein entsprechendes Programm erstellen und ablaufen lassen (→ Seite 71).
- Verwenden Sie den SmartShutdown-Assistenten (siehe unten), um das Programm zu erstellen und ablaufen zu lassen.

#### Erstellen eines Programms über den SmartShutdown-Assistenten

- 1. Öffnen Sie den Assistenten über SmartShutdown im Menü Batch.
- 2. Folgen Sie den Instruktionen auf den einzelnen Seiten des Assistenten. Klicken Sie auf die Schaltfläche **Hilfe**, wenn Sie weitere Informationen zu einer Seite benötigen.
- 3. Nach Beendigung des Assistenten
  - erstellt Chromeleon das Programm und speichert es unter der Zeitbasis ab, für die das Programm erstellt wurde.
  - öffnet Chromeleon das Dialogfenster Start Batch on.

Wählen Sie das Programm aus und starten Sie es mit Start.

Weitere Informationen zum SmartShutdown-Assistenten finden Sie in der *Chromeleon-Hilfe*.

# 5.9 Wartung und Wartungsintervalle

Der Autosampler ist aus hochwertigen Bauteilen und Materialien gefertigt und benötigt daher nur einen geringen Wartungsaufwand. Alle Oberflächen sind beständig gegen schwache Säuren, Basen und organische Lösungsmittel. Dennoch sollten Sie verschüttete oder verspritzte Flüssigkeiten sofort mit einem weichen, fusselfreien Tuch oder Papier aufsaugen (nicht trockenreiben). Eine längere Einwirkung kann Schäden verursachen.

Um die optimale Funktionsfähigkeit und maximale Verfügbarkeit Ihres Autosamplers sicherzustellen, führen Sie die folgenden Wartungsarbeiten in regelmäßigen Intervallen durch. Die Tabelle dient als Orientierungshilfe, welche Arbeiten Sie wann durchführen sollten. Wie häufig diese Arbeiten tatsächlich durchgeführt werden müssen, hängt von einer Vielzahl von Faktoren ab.

Häufigkeit	Was ist zu tun
Täglich	Prüfen Sie vor Betriebsbeginn die fluidischen Verbindungen auf Luftblasen. Entgasen Sie die Waschflüssigkeit (→ Seite 50).
	Prüfen Sie die Fluidik auf Undichtigkeiten.
	Prüfen Sie die fluidischen Verbindungen auf Salzablagerungen.
	Wenn Sie mit Pufferlösungen arbeiten, spülen Sie das System nach Abschluss der Arbeiten gründlich mit einer Flüssigkeit, die keine Puffer/Salze enthält.
Regelmäßig	Prüfen Sie die Schlauchverbindungen auf Beschädigungen (Knicke, Risse, Schnitte, Blockierung).
	Entfernen Sie die Probenhalter. Entfernen Sie auch die 5-Positionen Halter, falls nötig. Prüfen Sie den Drehteller auf Rückstande. Entfernen Sie Rückstände mit einem Tuch oder Papier. Diese Überprüfung ist speziell für Autosampler mit Proben- thermostatisierung wichtig, da ein sauberer Drehteller die Kühl- leistung verbessert.
	Prüfen Sie, ob an der Auslassöffnung des Injektionsventils Flüssigkeit austritt (→ Abb. 80, Seite 149). Abhilfemaßnahmen finden Sie auf Seite 149.
	Zur Ableitung von Flüssigkeit aus dem Innenraum sind an die beiden Abläufe rechts unterhalb des Autosamplers Schläuche angeschlossen (→ Seite 54). Prüfen Sie, dass die Schläuche nicht abgeknickt sind und an keiner Stelle höher als der Anschluss- Stutzen liegen. Entleeren Sie den Abfallbehälter, wenn erforderlich.
Häufigkeit	Was ist zu tun
------------	--
Jährlich	<ul> <li>Prüfen und tauschen Sie folgende Komponenten, falls erforderlich:</li> <li>Verbindungen am Injektionsventil (→ Seite 27)</li> <li>Verbindungen an der Spritze (→ Seite 53)</li> <li>Probennadel (→ Seite 125)</li> <li>Dichtung Injektionsventil (→ Seite 148)</li> </ul>
	<i>Empfehlung:</i> Lassen Sie den Kundendienst einmal pro Jahr eine vorbeugende Wartung durchführen.

Image: Image: Hinweis:Unter Chromeleon stehen Funktionen zur Lebensdauerprognose von<br/>Verschleißteilen sowie Diagnosetests für verschiedene Komponenten<br/>des Autosamplers zur Verfügung ( $\rightarrow$  Seiten 88 und 90).

# 6 Fehlersuche

## 6.1 Übersicht

Bei der Erkennung und Behebung von Fehlern, die beim Betrieb des Autosamplers oder eines UltiMate 3000-Systems auftreten können, helfen Ihnen:

## Status-LEDs

Die Status-LEDs (Light Emitting Diodes) auf der Gerätevorderseite lassen Sie auf den ersten Blick erkennen, ob der Autosampler eingeschaltet und mit Chromeleon verbunden ist. Darüber hinaus können Sie erkennen, ob der Autosampler korrekt arbeitet ( $\rightarrow$  Seite 20).

#### Meldungen

Wird während des Betriebs des Autosamplers ein Fehler erkannt, erscheint eine Meldung auf dem Gerätedisplay. Im Kapitel Meldungen auf dem Autosamplerdisplay finden Sie zu jeder Meldung Vorschläge für Abhilfemaßnahmen ( $\rightarrow$  Seite 104). Wird der Autosampler unter Chromeleon betrieben, wird der Fehler auch im Chromeleon Audit Trail protokolliert.

■ Hinweis: Hinweise zu Störungen, die beim Betrieb eines UltiMate 3000-Systems auftreten können, finden Sie im Kapitel Mögliche Störungen (→ Seite 112).

#### Diagnosetests

Wird der Autosampler unter Chromeleon betrieben, stehen Ihnen in Chromeleon eine Reihe von Testfunktionen zur Verfügung, mit denen Sie verschiedene Komponenten des Autosamplers auf ihre Funktionsfähigkeit hin testen können ( $\rightarrow$  Seite 89). Wurde ein Test nicht bestanden, finden Sie im Kapitel Diagnose-Meldungen Informationen zu möglichen Ursachen sowie Vorschläge für Abhilfemaßnahmen ( $\rightarrow$  Seite 110).

Wenn Sie einen Fehler mit Ihren Mitteln nicht beheben können, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.

## 6.2 Meldungen auf dem Autosamplerdisplay

Tritt während des Betriebs des Autosamplers ein Fehler auf, leuchtet die LED **Status** auf der Gerätevorderseite rot. Es erscheinen eine oder mehrere Meldungen auf dem Gerätedisplay. In der Navigationsleiste erscheinen die Funktionstasten **Prev**, **Next** und **Clear**.

Um	Wählen Sie
zur vorherigen Meldung zurück zu gelangen.	Prev
zur nächsten Meldung weiter zu gehen.	Next
die Meldung vom Gerätedisplay zu löschen.	Clear

Diese Tasten sind auch aktiv, wenn der Autosampler in Chromeleon verbunden ist.

Wenn der Autosampler in Chromeleon verbunden ist,

- wird der Fehler auch im Chromeleon Audit Trail protokolliert. Einige der genannten Meldungen sind für die verschiedenen Antriebe identisch. Im Chromeleon Audit Trail erscheint daher vor der eigentlichen Meldung der Antrieb, auf den sich die Meldung bezieht: Carousel (Drehteller), Syringe Drive (Spritzenantrieb), Needle Drive (Nadelantrieb) oder Horizontal Drive (horizontaler Antrieb). In der Tabelle haben diese Meldungen folgendes Format: (Antrieb): Meldung.
- können Sie Meldungen auf dem Gerätedisplay auch über den Chromeleon-Befehl ClearDisplayError löschen.

Die Tabelle listet Meldungen auf, die beim Betrieb des Autosamplers auftreten können, und nennt entsprechende Abhilfemaßnahmen. Zusätzlich können noch andere Meldungen erscheinen. Wenn Sie das Problem nicht beseitigen können, notieren Sie sich den genauen Wortlaut der Meldung. Wenden Sie sich an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.

Meldung	Abhilfe
24V power failure.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): End point not reached.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Error updating motor controller firmware.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Home position not found.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Home sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.

Meldung	Abhilfe
(Antrieb): Needle sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): No communication.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Not initialized.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Not ready.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Obstruction detected.	<ul> <li>Am Niederhalter befindet sich ein Hindernis. Entfernen Sie das Hindernis.</li> <li>Die Probennadel sticht auf den Rand des Probengefäßes, nicht in das Septum. Prüfen Sie, ob die Angaben auf der Registerkarte Segments / Pump Link in Chromeleon (→ Seite 41) mit der tatsächlichen Bestückung des Drehtellers übereinstimmt.</li> </ul>
(Antrieb): Stop condition at start.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
A self test is required.	Sie müssen einen Selbsttest durchführen. Der Autosampler funktioniert erst wieder, wenn der Selbsttest erfolgreich war. Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch. Wird während des Selbsttests ein Fehler gefunden, erscheint eine Meldung. Führen Sie die Abhilfemaßnahmen zu dieser Meldung durch.
Buffer loop is too full.	Die Pufferschleife enthält zu viel Flüssigkeit. Um die Pufferschleife zu leeren, geben Sie den <b>Buffer loop wash</b> im Menü <b>Control</b> $(\rightarrow$ Seite 78).
Carousel temperature sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Command input overflow.	Führen Sie ein Firmware- oder Treiberupdate durch.
Command queue overflow.	Das anwenderdefinierte Programm ist zu lang. Kürzen Sie das Programm.
Command syntax error.	Führen Sie ein Firmware- oder Treiberupdate durch.
Command was aborted.	Der Anwender hat den Befehl abgebrochen. Führen Sie den Befehl gegebenenfalls erneut aus.
Coolant temperature sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Cooling system overheated.	Vergewissern Sie sich, dass die Lüftungsschlitze des Autosamplers frei von Hindernissen sind. Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.

Meldung	Abhilfe
DVC position error.	Prüfen Sie, dass das Injektionsventil nicht blockiert ist. Die Überwurfmutter am Ventil ist gegebenenfalls zu fest angezogen. Wenn dies der Fall ist, kann das Ventil nicht mehr geschaltet werden. Lösen Sie die Überwurfmutter etwas. Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
	Vergewissern Sie sich, dass der Ventilkopf (Pod) korrekt installiert ist. Installieren Sie den Pod neu, falls erforderlich ( $\rightarrow$ Seite 150).
DVC (+ ergänzender Text) <i>oder</i> Error updating I2C controller firmware.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein und versuchen Sie es erneut. Installieren Sie die Firmware neu oder führen Sie gegebenenfalls ein Firmware-Update durch.
EEprom serialization error.	Führen Sie ein Firmware-Update durch.
Heat sink temperature sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Illegal command parameter.	Führen Sie ein Firmware- oder Treiberupdate durch.
Illegal syringe speed.	Die Spritzengeschwindigkeit ist nicht zulässig. Ändern Sie die Spritzengeschwindigkeit.
Illegal value.	Führen Sie ein Firmware- oder Treiberupdate durch.
Inject port missed.	Prüfen Sie, ob der Nadelsitz eingebaut ist. Installieren oder tauschen Sie den Nadelsitz, falls erforderlich ( $\rightarrow$ Seite 139).
	Prüfen Sie, ob die Probennadel eingebaut und intakt ist. Tauschen Sie die Probennadel, falls erforderlich ( $\rightarrow$ Seite 125).
	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Internal update. nn% finished. Do not interrupt. (wobei nn = Wert)	Es wird ein internes Update durchgeführt. Die Prozentzahl gibt den Grad der Ausführung an. Warten Sie, bis das interne Update beendet ist.
Invalid mix target.	Das Ziel für die Mischoperation ist nicht gültig. Geben Sie ein gültiges Ziel ein (Chromeleon-Befehle <b>UdpDraw &gt; From</b> und <b>UdpDispense &gt; To</b> ).
Leak detected.	Der Leaksensor des Autosamplers hat eine Undichtigkeit erkannt. Prüfen Sie, an welcher Stelle Flüssigkeit austritt, Beheben Sie die Ursache. Trocknen Sie den Leaksensor und die Auffangwanne (→ Seite 120).
Motor switching valve malfunction.	Prüfen Sie die Überwurfmutter am Injektionsventil. Wenn die Überwurfmutter zu fest angezogen ist, kann das Ventil nicht schalten. Lösen Sie die Überwurfmutter etwas. Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Motor switching valve timeout.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Needle home position not found.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.

Meldung	Abhilfe
Needle home sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Needle malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Property is protected.	Sie haben versucht, eine geschützte Variable zu ändern. Einige Variablen sind geschützt; der Anwender kann sie nicht ändern.
Property is read only.	Sie haben versucht, eine Lesevariable zu ändern. Einige Variablen sind Lesevariablen; der Anwender kann sie nicht ändern.
Reagent vial not found.	An der angegebenen Position befindet sich kein Reagensfläschchen. Stellen Sie ein Reagensfläschchen an die Position oder geben Sie eine andere Position an.
Remote control not enabled.	Die Verbindung zum Autosampler wurde in Chromeleon getrennt. Stellen Sie die Verbindung wieder her ("Connect").
	Führen Sie ein Firmware- oder Treiberupdate durch.
Sample height out of reach.	Bei dem installierten Probengefäß kann die Probennadel die ein- gestellte Sample Height nicht erreichen. Wählen Sie ein geeignetes Probengefäß oder ändern Sie die Einstellung für die Nadelhöhe $(\rightarrow$ Seite 91).
Syntax error.	Führen Sie ein Firmware- oder Treiberupdate durch
Syringe is not empty.	Zu Beginn der Injektion ist die Spritze nicht leer. Um die Spritze zu leeren, schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Syringe is too empty.	Es soll mehr Volumen aus der Spritze ausgestoßen werden, als zuvor aufgezogen wurde. Ändern Sie das Programm (auszustoßendes Volumen ≤ aufgezogenes Volumen).
Syringe is too full.	Die Spritze ist zu voll. Sie kann das definierte Ansaugvolumen nicht mehr aufnehmen. Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Syringe valve malfunction.	Prüfen Sie das Spritzenventil. Vergewissern Sie sich, dass die Anschlüsse am Ventil nicht zu fest angezogen sind.
Temperature difference too large. Probable coolant circuit malfunction.	Wenn der Drehteller offen ist, schließen Sie den Deckel. Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
The device is busy.	Das Gerät führt gerade ein anderes Kommando aus. Warten Sie, bis die Ausführung beendet ist. Führen Sie das Kommando erneut aus.
The volume is too large.	Es soll mehr Volumen aufgezogen werden, als die Spritze fassen kann. Ändern Sie das Volumen.
Unknown command.	Führen Sie ein Firmware- oder Treiberupdate durch
Unknown property.	Führen Sie ein Firmware- oder Treiberupdate durch.

Meldung	Abhilfe	
Unknown vial position.	Die angegebene Position existiert in der aktuellen Konfiguration nicht. Prüfen und korrigieren Sie die Segmenteinstellungen, falls erforderlich (→ Registerkarte Segments / Pump Link, Seite 41).	
USB interface is not ready.	Die USB-Verbindung konnte nicht hergestellt werden. Prüfen Sie die Verbindung. Wenn ein Kabel fehlerhaft ist, ersetzen Sie das Kabel.	
Vial not found.	An der angegebenen Position befindet sich kein Probengefäß. Stellen Sie ein Probengefäß an die angegebene Position oder geben Sie eine andere Position an.	
	Diese Meldung erscheint auch, wenn Ihre Probengefäße nicht die erforderliche Mindesthöhe haben ( $\rightarrow$ Seite 24). Verwenden Sie die Probengefäße, die auf Seite 24 genannt sind.	
	Es ist ein falscher Segmenttyp definiert. Prüfen und korrigieren Sie die Segmenteinstellungen, falls erforderlich ( $\rightarrow$ Registerkarte <b>Segments / Pump Link</b> , Seite 41).	
Vial pusher malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräte- rückseite aus und wieder ein oder führen Sie einen Selbsttest durch.	
Vial pusher obstructed.	Es befindet sich ein Hindernis am Niederhalter. Entfernen Sie das Hindernis.	
Wash port missed.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.	
Wrong tray type detected.	Die automatische Probenhaltererkennung hat festgestellt, dass die Angaben auf der Registerkarte <b>Segments</b> / <b>Pump Link</b> nicht mit der tatsächlichen Bestückung des Drehtellers übereinstimmen. Prüfen und korrigieren Sie die Angaben auf der Registerkarte <b>Segments</b> / <b>Pump</b> Link in Chromeleon (→ Seite 41).	
	Diese Meldung erscheint auch, wenn Ihre Probengefäße nicht die erforderliche Mindesthöhe (→ Seite 24) haben. Verwenden Sie die Probengefäße, die auf Seite 24 genannt sind.	

Wenn der Autosampler unter Chromeleon betrieben wird und keine Kommunikation zwischen dem Autosampler und Chromeleon aufgebaut werden kann, erscheinen Meldungen im Chromeleon Audit Trail.

Meldung	Abhilfe	
WPS-3000@USB-1610103 - Device not found on the USB.	Die USB-Verbindung zwischen dem Autosampler und dem Chromeleon-Server ist unterbrochen. Prüfen Sie die USB-Verbindung.	
	Die Stromversorgung des Autosamplers ist unterbrochen. Prüfen Sie die Verbindung des Autosamplers zum Stromnetz.	
Error opening WPS-3000 @USB-1610103 – The System cannot find the file specified	Die USB-Verbindung zwischen dem Autosampler und dem Chromeleon-Server ist unterbrochen. Prüfen Sie die USB-Verbindung. Die Stromversorgung des Autosamplers ist unterbrochen. Prüfen Sie die Verbindung des Autosamplers zum Stromnetz.	
Error issuing control request to WPS -3000@USB-1610103	Der Chromeleon-Server kann keine Verbindung zum Autosampler aufnehmen. Prüfen Sie die USB-Verbindung. Prüfen Sie die Verbindung des Autosamplers zum Stromnetz. Löschen Sie den in der Meldung genannten Autosampler aus der Konfiguration. Oder wählen in der Serverkonfiguration einen anderen Autosampler aus.	
Error reading from WPS-3000 @USB-1610103 Data error (cyclic redundancy check)	Prüfen Sie die USB Verbindung; die USB-Kabellänge darf 5 m zum nächsten Hub nicht überschreiten. Die maximale Gesamtkabellänge einschließlich der Hub-Verbindungen darf 30 m nicht überschreiten. Tauschen Sie fehlerhafte USB-Kabel aus. Tauschen einen fehlerhaften Hub aus.	
Error reading from WPS-3000 @USB-1610103	Die USB-Verbindung zwischen dem Autosampler und dem Chromeleon-Server ist unterbrochen. Prüfen Sie die USB-Verbindung. Die Stromversorgung des Autosamplers ist unterbrochen. Prüfen Sie die Verbindung des Autosamplers zum Stromnetz.	

## 6.3 Diagnose-Meldungen in Chromeleon

Wenn der Autosampler einen Diagnose-Test nicht besteht, führen Sie die genannten Abhilfemaßnahmen durch und wiederholen Sie den Test. Wird der Test auch dann nicht erfolgreich abgeschlossen, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.

#### **Drives and Sensors Test**

Wenn der Autosampler den Test nicht besteht, wiederholen Sie den Test. Schlägt der Test erneut fehl, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.

Testergebnis	Mögliche Ursache	Abhilfe
Needle seat leak too large.	Zu große Undichtigkeit im Injektionspfad (zwischen Injektionsventil, Probenschleife, Probennadel, Nadelsitz, Nadel- sitzkapillare und Injektionsventil).	<ul> <li>Prüfen Sie alle fluidischen Verbindungen im Injektionspfad auf Dichtigkeit.</li> <li>Prüfen Sie den Nadelsitz und die Probennadel.</li> <li>Tauschen Sie die den Nadelsitz, (→ Seite 125), die Nadelsitzkapillare (→ Seite 142) und/oder die Probennadel (→ Seite 139).</li> </ul>
System leak too large.	Zu hohe Undichtigkeit im System. Entweder ist die Pumpe undicht oder es besteht eine Undichtigkeit zwischen der Pumpe und dem Injektionsventil des Autosamplers.	<ul> <li>Prüfen Sie alle fluidischen Verbindungen zwischen der Pumpe und dem Injektionsventil im Autosampler auf Dichtigkeit.</li> <li>Testen Sie die Pumpe auf Dichtigkeit (→ Diagnosetests in Chromeleon oder <i>Bedienungsanleitung zur Pumpe</i>).</li> <li>Tauschen Sie das Rotor-Seal (oder Rotor-Seal und Stator Face Seal), den Ventilstator oder den Ventil-Pod aus (→ Kapitel 7.8.2 - 7.8.3, ab Seite 150).</li> </ul>

#### Leak Test fehlgeschlagen

Testergebnis	Mögliche Ursache	Abhilfe
The target pressure has not been reached, neither in INJECT nor in LOAD position.	Weder in der Position INJECT noch in der Position LOAD wurde der Zieldruck erreicht. Es gibt eine erhebliche Undichtigkeit im System.	<ul> <li>Prüfen Sie alle fluidischen Verbindungen auf Dichtigkeit.</li> <li>Prüfen Sie, ob an der Auslassöffnung des Injektionsventils Tröpfchen austreten (→ Abb. 80, Seite 149).</li> <li>Tauschen Sie das Rotor-Seal (oder Rotor-Seal und Stator Face Seal), den Ventilstator oder den Ventil-Pod aus (→ Kapitel 7.8.2 - 7.8.3, ab Seite 150).</li> <li>Testen Sie die Pumpe auf Dichtigkeit (→ Diagnosetests in Chromeleon</li> </ul>
		oder <i>Bedienungsanleitung zur</i> <i>Pumpe</i> ).
The target pressure has not been reached in the INJECT position.	In der Position INJECT wurde der eingestellte Druck nicht erreicht. Es gibt eine Undichtigkeit im Autosampler.	<ul> <li>Prüfen Sie alle fluidischen Verbindungen auf Dichtigkeit.</li> <li>Prüfen Sie speziell den Nadelsitz auf Anzeichen von Undichtigkeit.</li> <li>Prüfen Sie die fluidische Verbindung zwischen der Probennadel und der Probenschleife auf Undichtigkeiten.</li> <li>Prüfen Sie, ob an der Auslassöffnung des Injektionsventils Tröpfchen austreten (→ Abb. 80, Seite 149).</li> <li>Tauschen Sie das Rotor-Seal (oder Rotor-Seal und Stator Face Seal), den Ventilstator oder den Ventil-Pod aus (→ Kapitel 7.8.2 - 7.8.3, ab Seite 150).</li> </ul>

## 6.4 Mögliche Störungen

In der Tabelle finden Sie Hinweise zu Störungen, die beim Betrieb des UltiMate 3000-Systems auftreten können, deren mögliche Ursachen sowie entsprechende Abhilfemaßnahmen.

Weitere Hinweise und Abhilfemaßnahmen finden Sie in den Handbüchern zu den anderen Modulen eines UltiMate 3000-Systems.

Störung	Mögliche Ursache	Abhilfe
Keine Anzeige im Display.	Der Autosampler ist nicht am Stromnetz angeschlossen.	Schließen Sie das Netzkabel an.
	Der Netzschalter ist ausgeschaltet.	Schalten Sie den Netzschalter ein.
	Der Autosampler ist im Standby-Modus.	Drücken Sie die Taste <b>Standby</b> auf der Gerätevorderseite.
	Helligkeit und/oder Kontrast des Displays ist falsch eingestellt.	Stellen Sie die Helligkeit bzw. den Kontrast ein (→ Seite 86).
	Die Sicherung ist defekt.	Ersetzen Sie die Sicherungen $(\rightarrow \text{Seite 158}).$
	Die Ersatzsicherung brennt sofort durch.	Wenden Sie sich an den Kundendienst.
	Fehler in der Elektronik.	Wenden Sie sich an den Kundendienst.
Der Autosampler arbeitet nicht korrekt unter Chromeleon.	Es besteht keine Verbindung zwischen dem Autosampler und dem Chromeleon-Rechner.	Prüfen Sie das USB-Kabel und die Verbindung zum Chromeleon-Rechner.
	Die USB-Schnittstelle am Rechner ist nicht betriebsbereit	Prüfen Sie die USB- Schnittstelle am Rechner.
Kein Fluss	Im System tritt eine Undichtigkeit auf.	Beheben Sie die Ursache für die Undichtigkeit.
	Die verwendeten Eluenten sind alt oder verunreinigt oder die Qualität ist nicht ausreichend.	Tauschen Sie den Eluenten aus. Verwenden Sie Eluenten in HPLC-Qualität.
	Es gibt Verunreinigungen im System.	Spülen Sie das System mit einem geeigneten Lösungs- mittel.
	Weitere Ursachen finden Sie in der Bedienungsanleitung Ihrer Pumpe.	$\rightarrow$ Pumpenhandbuch
Im System herrscht ein hoher Gegendruck.	Kapillaren im System sind blockiert oder abgeknickt.	Prüfen Sie die Kapillaren im System schrittweise vom Detektor zur Pumpe. Tauschen Sie die Kapillaren, wenn erforderlich.

Störung	Mögliche Ursache	Abhilfe
Im System herrscht ein hoher Gegendruck <i>(Fortsetzung)</i>	Weitere Ursachen finden Sie in der Bedienungsanleitung Ihres Säulenthermostaten.	$\rightarrow$ TCC-Handbuch
Starke Basisliniendrift	Die Trennsäule ist verschmutzt.	Spülen oder ersetzen Sie die Säule.
	Das System ist nicht ausreichend äquilibriert.	Spülen Sie das System, bis Gleichgewicht erreicht ist. In der Regel ist dies nach 5-10 Säulenvolumina der Fall.
	Die verwendeten Eluenten sind verunreinigt oder nicht homogen.	Stellen Sie sicher, dass die Eluenten vor der Analyse vollständig durchmischt sind. Tauschen Sie den Eluenten aus. Prüfen Sie die Eluentenfilter. Achten Sie bei wässrigen Eluenten auf Verunreinigung durch Mikroorganismen.
	Die Umgebungsbedingungen sind instabil.	Stellen Sie eine gleichmäßige Temperatur und Luftfeuchtig- keit sicher. Vermeiden Sie Zugluft. Prüfen Sie am Detektor, dass die Abdeckungen der Lampen und Messzelle korrekt installiert sind und die Front- klappe geschlossen ist.
	Die mobile Phase wird im Kreislauf gefördert.	Fördern Sie die mobile Phase direkt in ein Abfallgefäß.
	Weitere Ursachen finden Sie in der Bedienungsanleitung Ihres Detektors.	$\rightarrow$ Detektorhandbuch
Starkes Rauschen, unregelmäßige Schwankungen der Basislinie	Der Eluent ist verunreinigt oder die Qualität ist nicht ausreichend.	Tauschen Sie den Eluenten aus. Verwenden Sie Eluenten in HPLC-Qualität.
	Weitere Ursachen finden Sie in der Bedienungsanleitung Ihres Detektors.	$\rightarrow$ Detektorhandbuch
Periodische Schwankungen der Basislinie, Pulsation	In der Pumpe treten Druck- schwankungen auf.	Entlüften und prüfen Sie die Pumpe (→ <i>Pumpenhandbuch</i> ).
	Es gibt Luftblasen im System.	Entlüften Sie das System $(\rightarrow Pumpenhandbuch)$ .
Peak-Tailing	Das Extrasäulenvolumen ist zu groß.	Verwenden Sie kurze Kapillaren mit geeignetem Innendurchmesser.

Störung	Mögliche Ursache	Abhilfe
Peak Tailing (Fortsetzung)	Die Kapillarverbindungen sind schlecht.	Tauschen Sie die Kapillaren aus. Verwenden Sie gegeben- enfalls Viper-Kapillaren.
Peakverbreiterung, hohe Totzeit	Es wird eine Kapillare mit einem zu großen Innendurch- messer verwendet.	Verwenden Sie eine Kapillare mit einem geeigneten Innen- durchmesser.
	Die Kapillaren sind blockiert oder die Kapillarverbindungen sind schlecht.	Tauschen Sie die Kapillaren aus. Verwenden Sie gegeben- enfalls Viper-Kapillaren.
	Der Eluentenfilter ist blockiert.	Prüfen Sie den Eluentenfilter auf Durchlässigkeit; tauschen Sie gegebenenfalls die Filter- fritten aus ( $\rightarrow$ <i>Pumpen-</i> <i>handbuch</i> ).
	Die Probenschleife ist verstopft.	Tauschen Sie die Proben- schleife aus (→ Seite 131).
	Das Proportionierventil der Pumpe ist defekt.	Wenden Sie sich an den Kundendienst.
	Die Trennsäule ist überladen oder verschmutzt.	Reinigen oder ersetzen Sie die Säule.
	Der Eluent hat sich verändert.	Verwenden Sie einen neuen Eluenten.
	Weitere Ursachen finden Sie in der Bedienungsanleitung Ihres Detektors.	$\rightarrow$ Detektorhandbuch
Reproduzierbar auftretende Störpeaks im Chromatogramm	Die Degaserkanäle sind verschmutzt.	Spülen Sie die Degaserkanäle (→ Handbuch zum Solvent Rack oder zur Pumpe).
	Die verwendeten Eluenten sind alt oder verunreinigt oder die Qualität ist nicht ausreichend.	Tauschen Sie den Eluenten aus. Verwenden Sie Eluenten in HPLC-Qualität.
	Es treten Verschmutzungen an anderen Stellen im System auf.	Spülen Sie das System mit einem geeigneten Lösungs- mittel.
Zusätzliche Peaks im Injektionspeak	Bei Gradienten ist die Äquili- brierzeit nach dem Spülschritt zu kurz.	Verlängern Sie die Äquili- brierzeit.
	Das Totvolumen ist zu hoch.	Beseitigen Sie das Totvolumen.
Einzelne breitere Störpeaks im Chromatogramm	Verspätete Elution aus einer vorausgegangenen Analyse.	Verlängern Sie die Laufzeit. Erhöhen Sie die Elutionsstärke des Gradienten (höherer organischer Anteil). Spülen Sie die Säule nach dem Probenlauf.

Störung	Mögliche Ursache	Abhilfe
Spikes	Es treten elektrische Störungen von anderen Geräten auf.	Isolieren Sie die Stromzufuhr von den anderen Geräten. Installieren Sie gegebenenfalls eine unterbrechungsfreie Stromversorgung (USV).
	Die Säulentemperatur liegt deutlich über dem Siedepunkt der mobilen Phase.	Installieren Sie einen Restriktor am Ausgang der Messzelle. Verwenden Sie einen Post- Column Cooler, wenn möglich $(\rightarrow TCC$ -Handbuch).
	Weitere Ursachen finden Sie in der Bedienungsanleitung Ihres Detektors.	$\rightarrow$ Detektorhandbuch
Negative Peaks	Lösungsmittel der Probe und mobile Phase unterscheiden sich in der Zusammensetzung.	Lösen Sie die Probe in der mobilen Phase.
	Die Absorption der gelösten Substanz ist geringer als die der mobilen Phase.	Messen Sie auf einer anderen Wellenlänge. Verwenden Sie eine mobile Phase mit geringerer Absorption.
Hohe Verschleppung	Die Nadelsitzkapillare ist nicht korrekt montiert. Dies kann zu einem erhöhten Totvolumen führen.	Prüfen Sie die Verbindung zwischen Nadelsitz und Nadelsitzkapillare. Wechseln Sie die Nadelsitzkapillare (→ Seite 142). Führt dies nicht zum Erfolg, müssen Sie auch den Nadelsitz tauschen (→ Seite 139).
	Die Probennadel ist verschlissen.	Tauschen Sie die Probennadel $(\rightarrow$ Seite 125).
Schlechte Reproduzierbarkeit bei großen Volumen	Im Probengefäß herrscht ein zu hoher Unterdruck.	Setzen Sie in Chromeleon den Parameter <b>PunctureOffset</b> . Der Drehteller wird um den einge- gebenen Wert gedreht, während sich die Probennadel im Septum befindet. Das Einstichloch wird entsprechend größer, so dass ein Druckausgleich stattfindet. Verwenden Sie keine Septen oder verwenden Sie geschlitzte Septen.

Störung	Mögliche Ursache	Abhilfe
Schlechte Präzision der Peakflächen	Der Autosampler saugt Luft an.	Es ist zu wenig Probe vor- handen, die Nadelhöhe ist nicht korrekt eingestellt (→ Seite 91) oder es gibt zu viele Replikate.
	Es gibt eine Luftblase in der Spritze.	Spülen Sie die Spritze (→ Seite 55).
	Es gibt eine Luftblase im Fluss- weg.	Führen Sie einen Wasch- vorgang durch (→ Seite 78). Die Waschflüssigkeit ist nicht entgast. Entgasen Sie die Waschflüssigkeit (→ Seite 50).
	Die Ansauggeschwindigkeit ist zu hoch.	Wählen Sie eine niedrigere Ansauggeschwindigkeit ( <b>DrawSpeed</b> ).
	Der Gasgehalt der Probe ist zu hoch oder gesättigt.	Wählen Sie eine niedrigere Ansauggeschwindigkeit ( <b>DrawSpeed</b> ). Entgasen Sie die Probe.
	Die Probennadel ist verstopft oder die Nadelspitze ist verformt.	Tauschen Sie die Probennadel aus ( $\rightarrow$ Seite 125).
	Der Autosampler, das Injektionsventil oder das Spritzenventil sind undicht.	Prüfen Sie den Autosampler und die Anschlüsse am Injektions- und Spritzenventil. Ziehen Sie undichte Verbin- dungen nach. Eventuell müssen das Injektionsventil und/oder Spritzenventil getauscht werden. Informationen zum Injektionsventil finden Sie im Kap. 7.8 (→ Seite 148). Wenn das Spritzenventil getauscht werden muss, wenden Sie sich an den Kundendienst.
	Es tritt Verschleppung auf.	Spülen Sie die Nadel mit einem geeigneten Lösungsmittel.
	Die Kapillarverbindungen sind nicht korrekt befestigt oder undicht.	Kontrollieren und ziehen Sie die Fittingverbindungen nach. Tauschen Sie den Nadelsitz (→ Seite 139).

Störung	Mögliche Ursache	Abhilfe
Schlechte Präzision der Peakflächen (Fortsetzung)	Es gibt Totvolumina in den Kapillarverbindungen.	Ersetzen Sie die Fitting- verbindungen. Achten Sie auf korrekte Installation der Verbindungen. Verwenden Sie gegebenenfalls Viper- Kapillaren.
	Die Kolbendichtringe sind undicht.	Tauschen Sie die Dichtungen aus ( $\rightarrow$ <i>Pumpenhandbuch</i> ).
	Es befindet sich Luft im Arbeitskolben.	Entlüften und prüfen Sie die Pumpe (→ <i>Pumpenhandbuch</i> ).
	Die Pumpe pulsiert.	Verwenden Sie entgastes Lösungsmittel.
	Der Gradient ist nicht reproduzierbar.	Ändern Sie den Gradienten. Prüfen Sie die Pumpenfunktion und Entgasung. Prüfen Sie die Ansaugfritten auf Verstopfung. Tauschen Sie die Fritten.
	Die Probe ist instabil und zerfällt.	Verwenden Sie eine neue Probe oder ändern Sie die Bedin- gungen. Kühlen Sie die Probe im Autosampler, wenn möglich.
	Basislinienschwankungen	Siehe Abhilfen, die bei den Störungen zur Basislinie beschriebenen sind. (weiter oben in der Tabelle).
	Die Umgebungsbedingungen sind instabil.	Stellen Sie eine gleichmäßige Temperatur und Luft- feuchtigkeit sicher. Vermeiden Sie Zugluft. Verwenden Sie gegebenenfalls einen Säulen- thermostaten.
	Es treten Verschmutzungen im System auf.	Spülen Sie das System mit einem geeigneten Lösungsmittel.

# 7 Service

## 7.1 Allgemeine Hinweise und Sicherheitsmaßnahmen

Die nachfolgenden Abschnitte beschreiben die Service- und Reparaturarbeiten, die Sie als Anwender ausführen dürfen. Weitergehende Reparaturarbeiten dürfen nur vom Thermo Fisher Scientific-Kundendienst ausgeführt werden.

Warnung: Die fluidischen Komponenten des Gerätes können mit gesundheitsschädlichen Lösungsmitteln gefüllt sein. Ebenso können gesundheitsschädliche Substanzen an den medienberührten Teilen anhaften. Tragen Sie eine geeignete Schutzausrüstung. Spülen Sie die fluidischen Komponenten mit einem geeigneten Lösungsmittel von gesundheitsschädlichen Substanzen frei.

> Informationen zum richtigen Umgang mit konkreten Substanzen und Empfehlungen für konkrete Gefahrensituationen entnehmen Sie bitte dem Sicherheitsdatenblatt der Substanzen, mit denen Sie umgehen. Beachten Sie die Richtlinien der Guten Laborpraxis (GLP).

Bevor Sie Service- und Reparaturarbeiten ausführen, beachten Sie die folgenden Hinweise:

- Beachten Sie bei allen Servicearbeiten die in dieser Anleitung angegebenen Sicherheitshinweise.
- Vor Arbeiten an der Fluidik müssen Sie die Spritze und/oder Nadel in die geeignete Position fahren. Um Beschädigungen der Spritze und des Autosamplers zu vermeiden, legen Sie keine Gegenstände unter den Spritzenstößel.
- *Empfehlung* Testen Sie den Autosampler auf Dichtigkeit, nachdem Sie die Wartungsarbeiten an der Fluidik vorgenommen haben. Führen Sie in Chromeleon den **Leak Test** aus.
- Verwenden Sie ausschließlich die von Thermo Fisher Scientific für das Gerät autorisierten Original-Ersatzteile.
- Falls ein Gerät zur Reparatur zurückgeschickt werden muss, wenden Sie sich zunächst an den Thermo Fisher Scientific Kundendienst für Dionex HPLC-Produkte. Für die Rücksendung ist eine RMA- (Return Material Authorization) Nummer erforderlich. Der Transport darf nur in der Originalverpackung unter Beachtung der Verpackungsvorschrift erfolgen. Erfolgt die Einsendung nicht in der Originalverpackung, entfällt die Gerätegarantie.

Ist die Originalverpackung nicht mehr verfügbar, können Sie geeignete Geräteverpackungen über die Thermo Fisher Scientific-Vertriebsorganisation für Dionex HPLC-Produkte bestellen. Die Verpackungsvorschrift ist im Ordner "Installation and Qualification Documents for Chromatography Instruments" enthalten. Sie sind auch auf Anfrage erhältlich.

Hinweise zur Außerbetriebnahme des Autosamplers finden Sie auf Seite 98.

## 7.2 Beseitigen von Undichtigkeiten

Der Leaksensor spricht an, wenn er Flüssigkeit ausgesetzt ist. Beseitigen Sie die Ursache für die Undichtigkeit und trocknen Sie den Leaksensor:

- 1. Schalten Sie den Autosampler aus.
- 2. Prüfen Sie den Autosampler auf eventuelle Undichtigkeit. Ziehen Sie undichte Schlauch- und Kapillarverbindungen nach oder tauschen Sie diese aus.
- 3. Saugen Sie mit einem Tuch die Flüssigkeit auf, die sich am unteren Ende des Leaksensors in der Auffangwanne gesammelt hat.

**Vorsicht:** Seien Sie vorsichtig. Verbiegen oder beschädigen Sie die Sensoren nicht.



Abb. 37: Trocknen des Leaksensors

- 4. Lassen Sie dem Sensor einige Minuten Zeit, sich auf die Umgebungstemperatur einzustellen
- 5. Schalten Sie den Autosampler ein.
- 6. Wird nach dem Einschalten des Autosamplers kein Fehler gemeldet, können Sie den Betrieb wieder aufnehmen.
- **i Hinweis:** Die LED **Status** auf der Gerätevorderseite bleibt rot, bis der Sensor trocken ist. Wenn eine Meldung auf dem Gerätedisplay erschienen ist, lässt sich diese über die Taste **Clear** auf der Navigationsleiste löschen  $(\rightarrow$  Seite 77).

## 7.3 Wechseln der Spritze

Beschreibung	BestNr.
WPS-3000SL Analytical, WPS-3000TBSL Analytical, WPS-3000RS, WPS-3000TBRS:	
100-µL-Spritze	6822.0002
WPS-3000TXRS: 25-µL-Spritze	6822.0001
WPS-3000SL Semiprep: 2500-µL-Spritze	6822.0006
Informationen zu optional erhältlichen Spritzen finden Sie im Kapitel 10.2 ( $\rightarrow$ Seite 178).	

- Bewegen Sie die Spritze nach unten.
   Rufen Sie am Gerät das Menü Maintenance auf und wählen Sie Syringe und Down oder—geben Sie in Chromeleon das Kommando InitiateChangeSyringe.
- 2. Klappen Sie den Frontdeckel nach oben auf.
- Lösen Sie die Halteschraube f
  ür den Spritzenst
  ö
  ßel mit einem Maulschl
  üssel (Gr
  ö
  ße 8). Entfernen Sie die Schraube per Hand. Heben Sie die Schraube f
  ür den Einbau der neuen Spritze auf.



Abb. 38: Lösen der Spritze am Spritzenstößel

4. Schrauben Sie die Spritze am Spritzenventil ab.



Abb. 39: Lösen der Spritze am Spritzenventil

5. Vergewissern Sie sich, dass beim Ausbau der alten Spritze die Dichtung im Ventil geblieben ist. Schauen Sie zum Beispiel einem Spiegel in den unteren Anschlussport des Ventils.

Falls die Dichtung zusammen mit der alten Spritze entfernt wurde, installieren Sie eine neue Dichtung, bevor Sie die neue Spritze anschließen. Die Dichtung ist im Lieferumfang der Ersatzspritze enthalten; sie ist auch als separates Ersatzteil erhältlich (Best.-Nr. 6822.0009 für 5 Dichtungen).

**Vorsicht:** Im Spritzenventil darf sich nur eine Dichtung befinden. Wenn Sie zwei oder mehr Dichtungen installieren, wird das Spritzenventil beschädigt.

- 6. Entfernen Sie die Luft aus der neuen Spritze. Füllen Sie die neue Spritze mit einer Plastikspritze oder gehen Sie wie unten beschrieben vor:
  - a) Füllen Sie ein Becherglas mit der Waschflüssigkeit oder einer Mischung aus Isopropanol und Wasser (50:50).
  - b) Geben Sie die Spritze mit dem Spritzeneingang nach unten in das Becherglas.
  - c) Bewegen Sie die Spritze mit eingefahrenem Stößel mehrfach schnell hin und her. Damit werden Luftblasen entfernt, die sich unter Umständen am Spritzeneingang befinden.
  - d) Ziehen Sie die Spritze mehrfach hintereinander zügig auf und stoßen Sie den Inhalt schnell aus, bis die Spritze luftfrei ist.
- 7. Ziehen Sie die Spritze komplett auf.
- 8. Richten Sie die Spritze unter dem Ventil aus (das Gewinde zeigt nach oben).
- 9. Halten Sie die Spritze senkrecht und stoßen Sie etwas Flüssigkeit aus, damit sich am Spritzeneingang keine Luft mehr befindet. Schrauben Sie die Spritze am Spritzenventil fest.



Flüssigkeitsfilm

Abb. 40: Flüssigkeitsfilm am Spritzeneingang

- 10. Richten Sie den Spritzenstößel aus. Schrauben Sie die Befestigungsschraube mit der Hand ein. Ziehen Sie die Schraube mit dem Maulschlüssel etwas fester.
- 11. Klappen Sie den Frontdeckel nach unten.

12. Bewegen Sie die Spritze nach oben.

Rufen Sie am Gerät das Menü Maintenance auf und wählen Sie Syringe und Upoder —geben Sie in Chromeleon das Kommando TerminateChangeSyringe.

**Vorsicht:** Vergewissern Sie sich, dass die Spritze korrekt montiert ist, ehe Sie die Spritze nach oben fahren. Andernfalls kann die Spritze beschädigt werden.

- Spülen Sie die neue Spritze mehrmals (→ Kapitel 4.6, Seite 55).
   Luft in der Spritze kann die Messergebnisse stark beeinflussen. Bevor Sie eine Analyse starten, muss die Spritze luftfrei sei. Wenn in der Spritze noch Luftblasen zu sehen sind, wiederholen Sie den Spülvorgang.
- 14. Wenn die neue Spritze ein anderes Volumen hat als die Spritze, die Sie ausgebaut haben:
  - ♦ Prüfen Sie in der Konfiguration des Autosamplers auf der Seite Options, dass die Spritzengröße korrekt angegeben ist (→ Seite 43). Ändern Sie die Einstellung, falls erforderlich. Übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Klicken Sie auf der Seite General auf die Schaltfläche Send configuration to module (→ Seite 45).
  - ◆ Prüfen Sie, dass passende Puffer- und Probenschleifen installiert sind (→ Seite 17). Das Volumen der Probenschleife sollte mindestens dem Volumen der Spritze entsprechen.
- 15. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl **SyringeChanged**.

## 7.4 Wechseln der Pufferschleife

Beschreibung	BestNr.
WPS-3000SL Analytical, WPS-3000TBSL Analytical, WPS-3000RS, WPS-3000TXRS: Pufferschleife, 100 μL, SST, Viper	6820.2466
WPS-3000TBRS: Pufferschleife, 100 μL, MP35N,Viper	6841.2466
WPS-3000SL Semiprep: Pufferschleife, >250 μL, SST, Viper	6820.2468

- 1. Klappen Sie den Frontdeckel nach oben auf.
- 2. Die Pufferschleife ist mit einem Halteclip am Injektionsventilhalter befestigt. Öffnen Sie den Clip und entfernen Sie die Pufferschleife.
- 3. Schrauben Sie die Pufferschleife am Spritzenventil ab.



Schrauben Sie die Pufferschleife vom Spritzenventil ab.

Abb. 41: Pufferschleife am Spritzenventil

- 4. Entfernen Sie die Pufferschleife am Injektionsventil.
- 5. Befestigen Sie die neue Pufferschleife in der Halterung am Injektionsventilhalter.
- 6. Schrauben Sie die neue Pufferschleife am Injektionsventil und am Spritzenventil fest.
- 7. Klappen Sie den Frontdeckel nach unten.
- 8. Führen Sie einen Waschzyklus für die Pufferschleife durch.

Rufen Sie am Gerät das Menü **Control** auf und wählen Sie **Buffer loop wash**—*or*— geben Sie in Chromeleon das Kommando **WashBufferLoop**.

Das Waschvolumen sollte mindestens das doppelte Volumen der Pufferschleife sein. Ziehen Sie die Spritze so oft auf wie erforderlich.

- 9. Wenn die neue Pufferschleife ein anderes Volumen hat als die Pufferschleife, die Sie ausgebaut haben, vergewissern Sie sich, dass auch eine geeignete Probenschleife und Spritze installiert ist (→ Seite 17).
- 10. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

## 7.5 Probennadel

## 7.5.1 Wechseln der Probennadel

Die Schritte zum Wechseln der Probennadel hängen vom Autosamplermodell ab:

- WPS-3000SL Analytical, WPS-3000RS, WPS-3000TBRS, WPS-3000TXRS Folgen Sie den Schritten im Kapitel 7.5.1.1.
- WPS-3000TBSL Analytical und WPS-3000SL Semiprep Folgen Sie den Schritten im Kapitel 7.5.1.2 ( $\rightarrow$  Seite 128).

**Warnung:** Die Probennadel ist sehr spitz und gegebenenfalls kontaminiert. Seien Sie vorsichtig, um sich nicht zu verletzen. Spülen Sie die Probennadel von gesundheitsschädlichen Substanzen frei.

#### 7.5.1.1 WPS-3000SL Analytical, WPS-3000RS, WPS-3000TBRS, WPS-3000TXRS

Diese Kapitel gilt nicht für den WPS-3000TBSL Analytical. Für diesen Autosampler, siehe Kapitel 7.5.1.2 ( $\rightarrow$  Seite 128).

Beschreibung	BestNr.
WPS-3000SL Analytical, WPS-3000RS, WPS-3000TXRS: Probennadel, SST, Viper-kompatibel	6820.2432
WPS-3000TBRS Probennadel, Keramik, Viper-kompatibel	6841.2420

1. Bewegen Sie die Probennadel in die Position für den Nadelwechsel (Probennadel oben und Nadelarm ganz links).

Rufen Sie am Gerät das Menü Maintenance auf und wählen Sie Needle Position und Maintain— oder — geben Sie in Chromeleon das Kommando InitiateChangeNeedle.

2. Klappen Sie den Frontdeckel nach oben auf.

3. Eine beige Schraube fixiert die Einheit aus Probennadel und Probenschleife im Nadelhalter. Lösen Sie Schraube. Ziehen Sie die Einheit vorsichtig nach oben aus dem Nadelhalter heraus. Achten Sie darauf, die Probenschleife nicht zu verbiegen.

Beige Sicherungsschraube Probennadel



Abb. 42: Entfernen der Probennadel aus dem Nadelhalter

 Lösen Sie die Verbindung zwischen Probenschleife und Probennadel.
 Halten Sie die Probennadel mit dem Maulschlüssel fest (Größe 5.5 mm; im Zubehör enthalten). Lösen Sie die Verbindung per Hand mit der Rändelschraube.



Abb. 43: Lösen der Verbindung Probenschleife - Probennadel

Behalten Sie die Sicherungsschraube für den Einbau der neuen Probennadel.

- 5. Verbinden Sie die neue Probennadel und die Probenschleife. Achten Sie darauf, die Komponenten nicht zu verbiegen.
  - a) Vergewissern Sie sich, dass sich die beige Sicherungsschraube auf der Probenschleife befindet. Wenn nicht, schieben Sie die Schraube wieder auf die Schleife.



Abb. 44: Probenschleife mit Sicherungsschraube

- b) Schieben Sie die Probenschleife bis zum Anschlag in die Probennadel hinein.
- c) Halten Sie die Probennadel mit dem Maulschlüssel. Ziehen Sie die Verbindung per Hand mit der schwarzen Rändelschraube fest.
- d) Setzen Sie die Einheit in den Nadelhalter ein. Fixieren Sie die Einheit mit der Sicherungsschraube Probennadel.

- 6. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test** *oder* geben Sie in Chromeleon das Kommando **SelfTest**.
- 7. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl NeedleChanged.
- Empfehlung Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch (→ Seite 90).

**i Hinweis:** We chseln Sie nach der Probennadel auch der Nadelsitz ( $\rightarrow$  Seite 139).

## 7.5.1.2 WPS-3000TBSL Analytical und WPS-3000SL Semiprep

Beschreibung	BestNr.
WPS-3000TBSL Analytical: Probennadel, SST	6820.2403
WPS-3000SL Semiprep: Probennadel, SST	6820.2419

1. Bewegen Sie die Probennadel in die Position für den Nadelwechsel (Probennadel oben und Nadelarm ganz links).

Rufen Sie am Gerät das Menü **Maintenance** auf und wählen Sie **Needle Position** und **Maintain**— *oder*—geben Sie in Chromeleon das Kommando **InitiateChangeNeedle**.

- 2. Klappen Sie den Frontdeckel nach oben auf.
- 3. Eine beige Schraube fixiert die Einheit aus Probennadel und Probenschleife im Nadelhalter. Lösen Sie Schraube. Ziehen Sie die Einheit vorsichtig nach oben aus dem Nadelhalter heraus. Achten Sie darauf, die Probenschleife nicht zu verbiegen.



Abb. 45: Entfernen der Probennadel aus dem Nadelhalter (links WPS-3000SL Semiprep; rechts WPS-3000TBSL Analytical)

4. Lösen Sie die Verbindung zwischen Probenschleife und Probennadel mit den beiden Maulschlüsseln aus dem Zubehör.





Abb. 46: Lösen der Verbindung Probenschleife - Probennadel (links: WPS-3000SL Semiprep; rechts: WPS-3000TBSL Analytical)

- 5. Bauen Sie die neue Probennadel in umgekehrter Reihenfolge ein. Zur leichteren Installation sollten Sie die Verbindung zwischen Probennadel und Probenschleife zu zweit herstellen. Schieben Sie die Probenschleife bis zum Anschlag in die Probennadel hinein. Halten Sie beide Teile in dieser Position fest, während Sie die Verbindung mit den Maulschlüsseln festziehen.
- 6. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder*—geben Sie in Chromeleon das Kommando **SelfTest**.
- 7. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl NeedleChanged.
- 8. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

**i** Hinweis: Wechseln Sie nach der Probennadel auch der Nadelsitz ( $\rightarrow$  Seite 139).

## 7.5.2 Waschen der Probennadel

Folgende Optionen stehen zum Waschen der Probennadel zur Verfügung:

- Spülen der Probennadel von außen Dabei wird das Äußere der Probennadel im Waschport mit Waschflüssigkeit gespült. Nach dem Waschen wird die Probennadel über die Membranpumpe mit Luft getrocknet.
- Spülen des Inneren der Probennadel Um Luft oder irrtümlich aufgezogene Probe aus der Probennadel zu entfernen, können Sie das Innere der Probennadel mit Waschflüssigkeit spülen.

## 7.5.2.1 Waschen der Probennadel von außen

#### Über Chromeleon

- 1. Öffnen Sie das Dialogfenster Commands.
- 2. Klicken Sie unter **Sampler** auf **WashVolume**. Geben Sie Volumen ein, mit dem der Waschport gespült werden soll.
- 3. Legen Sie unter **WashSpeed** die Geschwindigkeit der Spritze für den Waschvorgang fest.
- 4. Starten Sie den Waschvorgang mit dem Kommando Wash.

#### Vom Gerätedisplay aus

- Wählen Sie die Funktionstaste **Wash**. Dabei wird das Volumen verwendet, das Sie unter **Wash** im Menü **Control** festgelegt haben.
- Rufen Sie das Menü **Control** auf und wählen Sie **Wash**. Legen Sie das Waschvolumen fest und führen Sie den Waschzyklus durch.

#### 7.5.2.2 Waschen der Probennadel von innen

- Öffnen Sie in Chromeleon das Dialogfenster **Commands** und geben Sie den Befehl **WashSampleLoop**.
- Rufen Sie am Gerätedisplay das Menü Control auf und wählen Sie Inject Loop Wash.
- **i** Hinweis: Wenn Sie den Befehl WashSampleLoop oder Inject Loop Wash gegeben haben, wird bei der folgenden Injektion (Selbsttest oder InjectValveToInject-Befehl) Reinigungslösung injiziert. Dies kann die Äquilibrierung einer angeschlossenen Säule stören.

## 7.6 Wechseln der Probenschleife

Die Schritte zum Wechseln der Probenschleife hängen vom Autosamplermodell ab.

- WPS-3000SL Analytical, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS: Folgen Sie den Schritten im Kapitel 7.6.1.
- WPS-3000TBSL Analytical: Folgen Sie den Schritten im Kapitel 7.6.2 (→ Seite 134).
- WPS-3000SL Semiprep: Folgen Sie den Schritten im Kapitel 7.6.3 (→ Seite 136).

## 7.6.1 WPS-3000SL Analytical, WPS-3000RS/TBRS/TXRS

Dieses Kapitel gilt nicht für den WPS-3000TBSL Analytical. Für diesen Autosampler, siehe Kapitel 7.6.2 ( $\rightarrow$  Seite 134).

Beschreibung	BestNr.
<i>WPS-3000SL Analytical und WPS-3000RS</i> : Probenschleife, analytisch, 100 $\mu$ L (V= 130 $\mu$ L) <sup>1</sup> , SST, Viper	6820.2451
<i>WPS-3000TBRS</i> : Probenschleife, Mikro, 25 $\mu$ L (V= 40 $\mu$ L) <sup>1</sup> , MP35N, Viper	6841.2452
<i>WPS-3000TXRS</i> : Probenschleife, Mikro, 25 $\mu$ L (V= 40 $\mu$ L) <sup>1</sup> , SST, Viper	6820.2452
Informationen zu optional erhältlichen Probenschleifen finden Sie im Kapitel 10.2 (→ Seite 178).	

1 Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen.

- Bewegen Sie die Probennadel in die geeignete Position f
  ür den Wechsel der Probenschleife (Probennadel oben und Nadelarm ganz links). Rufen Sie am Ger
  ät das Men
  ü Maintenance auf und w
  ählen Sie Needle Position und Maintain— oder — geben Sie in Chromeleon das Kommando InitiateChangeNeedle.
- 2. Klappen Sie den Frontdeckel nach oben auf.

Warnung: Die Probenschleife ist mit der Probennadel verbunden. Die Probennadel ist sehr spitz und gegebenenfalls kontaminiert. Seien Sie vorsichtig, um sich nicht zu verletzen. Spülen Sie die Probennadel von gesundheitsschädlichen Substanzen frei.

3. Eine beige Schraube fixiert die Einheit aus Probennadel und Probenschleife im Nadelhalter. Lösen Sie Schraube. Ziehen Sie die Einheit vorsichtig nach oben aus dem Nadelhalter heraus. Achten Sie darauf, die Probenschleife nicht zu verbiegen.

Beige Sicherungsschraube (Probennadel)



Abb. 47: Entfernen der Probennadel aus dem Nadelhalter

4. Lösen Sie die Verbindung zwischen Probenschleife und Probennadel.

Halten Sie die Probennadel mit dem Maulschlüssel fest (Größe 5.5 mm, im Zubehör enthalten). Lösen Sie die Verbindung per Hand mit der schwarzen Rändelschraube.



Abb. 48: Lösen der Verbindung Probenschleife - Probennadel

Behalten Sie die Sicherungsschraube für den Einbau der neuen Probenschleife.

5. Die Probenschleife ist in einem Halteclip am Injektionsventilhalter befestigt. Entfernen Sie die Probenschleife mit dem Clip. (Die neue Probenschleife wird mit Clip und Befestigungsschraube geliefert.)

> Entfernen Sie die Probenschleife mit dem Clip.



Abb. 49: Halterung am Injektionsventil

- 6. Entfernen Sie die Probenschleife am Injektionsventil.
- 7. Verbinden Sie die Probennadel und die neue Probenschleife. Achten Sie darauf, die Komponenten nicht zu verbiegen.
  - a) Schieben Sie die beige Sicherungsschraube Probennadel auf die neue Probenschleife.



Abb. 50: Probenschleife mit Sicherungsschraube

- b) Schieben Sie die Probenschleife bis zum Anschlag in die Probennadel hinein.
- c) Halten Sie die Probennadel mit dem Maulschlüssel. Ziehen Sie die Verbindung per Hand und mit der schwarzen Rändelschraube fest.
- 8. Verbinden Sie die Probenschleife mit dem Injektionsventil.
- 9. Befestigen Sie die Probenschleife am Injektionsventilhalter ( $\rightarrow$  Abb. 49).
- 10. Setzen Sie die Einheit aus Probennadel und Probenschleife in den Nadelhalter ein. Fixieren Sie die Einheit mit der beigen Sicherungsschraube Probennadel.
- 11. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder* geben Sie in Chromeleon das Kommando **SelfTest**.
- 12. Wenn die neue Probenschleife ein anderes Volumen hat als die Probenschleife, die Sie ausgebaut haben:
  - ◆ Prüfen Sie in der Konfiguration des Autosamplers auf der Seite Options, dass s das Volumen der Probenschleife korrekt angegeben ist (→ Seite 43). Ändern Sie die Einstellung, falls erforderlich. Übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Klicken Sie auf der Seite General auf die Schaltfläche Send configuration to module (→ Seite 45).
  - ♦ Prüfen Sie, dass eine geeignete Pufferschleife und Spritze installiert ist (→ Seite 17). Das Volumen der Probenschleife sollte mindestens dem Volumen der Spritze entsprechen.
- 13. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

## 7.6.2 WPS-3000TBSL Analytical

Beschreibung	BestNr.
Probenschleife, analytisch, 100 $\mu$ L (V= 150 $\mu$ L) <sup>1</sup> , PEEK	6820.2431

1 Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen.

- 1. Bewegen Sie die Probennadel in die Position für den Nadelwechsel (Probennadel oben und Nadelarm ganz links). Rufen Sie am Gerät das Menü **Maintenance** auf und wählen Sie **Needle Position** und **Maintain** *oder* geben Sie in Chromeleon das Kommando **InitiateChangeNeedle**.
- 2. Klappen Sie den Frontdeckel nach oben auf.
- 3. Entfernen Sie die Probenschleife aus der Halterung am Injektionsventilhalter.



Abb. 51: Halterung am Injektionsventil

- 4. Entfernen Sie die Probenschleife am Injektionsventil.
- 5. Eine beige Schraube fixiert die Einheit aus Probennadel und Probenschleife im Nadelhalter. Lösen Sie Schraube. Ziehen Sie die Einheit vorsichtig nach oben aus dem Nadelhalter heraus.



Abb. 52: Entfernen der Probennadel aus dem Nadelhalter

6. Lösen Sie die Verbindung zwischen Probenschleife und Probennadel mit den beiden Maulschlüsseln aus dem Zubehör.



Abb. 53: Lösen der Verbindung Probenschleife - Probennadel

- Bauen Sie die neue Probennadel in umgekehrter Reihenfolge ein. Beachten Sie die folgenden Hinweise:
  - Zur leichteren Installation sollten Sie die Verbindung zwischen Probennadel und Probenschleife zu zweit herstellen. Schieben Sie die Probenschleife bis zum Anschlag in die Probennadel hinein. Halten Sie beide Teile in dieser Position fest, während Sie die Verbindung mit den Maulschlüsseln festziehen
  - ◆ Die grüne Hülse der Probenschleife muss in der Halterung am Injektionsventilhalter liegen (→ Abb. 54). Nur dann ist die Probenschleife korrekt installiert.



Abb. 54: Probenschleife, eingebaut

- 8. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder* geben Sie in Chromeleon das Kommando **SelfTest**.
- 9. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl NeedleChanged.
- 10. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

## 7.6.3 WPS-3000SL Semiprep

Beschreibung	BestNr.
Probenschleife, semiprep, 2500 $\mu$ L (V = 3300 $\mu$ L) <sup>1</sup> , SST	6820.2416
Informationen zu optional erhältlichen Probenschleifen finden Sie im Kap ( $\rightarrow$ Seite 178).	pitel 10.2

1 Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen.

 Bewegen Sie die Probennadel in die geeignete Position f
ür den Wechsel der Probenschleife (Probennadel oben und Nadelarm ganz links).
 Rufen Sie am Ger
ät das Men
ü Maintenance auf und w
ählen Sie Needle Position und Maintain— oder —geben Sie in Chromeleon das Kommando InitiateChangeNeedle.

- 2. Klappen Sie den Frontdeckel nach oben auf.
- 3. Lösen Sie die beiden Schrauben auf dem Halteblech für die Probenschleife mit circa 2 Umdrehungen (Schrauben nicht entfernen). Ziehen Sie die Probenschleife nach hinten aus dem Halteblech heraus.



Ziehen Sie die Probenschleife nach hinten heraus.

Lösen Sie diese beiden Schrauben.

Abb. 55: Halteblech Probenschleife

4. Entfernen Sie die Probenschleife aus der Halterung am Injektionsventilhalter.



Entfernen Sie die Probenschleife aus dieser Halterung.

Abb. 56: Halterung am Injektionsventil
5. Die Probenschleife wird in Halterungen an der inneren Rückwand entlang geführt. Nehmen Sie die Probenschleife nach oben aus diesen Halterungen heraus.



Abb. 57: Haltung für Probenschleife

- 6. Entfernen Sie die Probenschleife am Injektionsventil.
- 7. Eine Fittingschraube fixiert die Einheit aus Probennadel und Probenschleife im Nadelhalter. Lösen Sie Schraube. Ziehen Sie die Einheit vorsichtig nach oben aus dem Nadelhalter heraus.



Abb. 58: Entfernen der Probennadel aus dem Nadelhalter

8. Lösen Sie die Verbindung zwischen Probenschleife und Probennadel mit den beiden Maulschlüsseln aus dem Zubehör.



Abb. 59: Lösen der Verbindung Probenschleife - Probennadel

- 9. Bauen Sie die neue Probenschleife in umgekehrter Reihenfolge ein. Beachten Sie die folgenden Hinweise:
  - An den beiden Enden der Probenschleife befinden sich unterschiedliche Ferrulen:





Diese Ferrule muss sich an dem Ende befinden, dass mit der Probennadel verbunden wird.

Abb. 60: Unterschiedliche Ferrulen bei einer semipräparativen Probenschleife

• Schieben Sie das Fitting auf die Probenschleife (Nadelende), ehe Sie die Probenschleife mit der Probennadel verbinden.



Schieben Sie das Fitting auf das Ende der Probenschleife, dass mit der Probennadel verbunden wird.

Abb. 61: Fitting für das Nadelende der Probenschleife

• Zur leichteren Installation sollten Sie die Verbindung zwischen Probennadel und Probenschleife zu zweit herstellen. Die Probennadel ist sehr spitz. Gehen Sie vorsichtig vor, um sich nicht zu verletzten

Schieben Sie die Probenschleife bis zum Anschlag in die Probennadel hinein. Halten Sie beide Teile in dieser Position fest, während Sie die Verbindung mit den Maulschlüsseln festziehen.

- 10. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück .Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder* geben Sie in Chromeleon das Kommando **SelfTest**.
- 11. Wenn die neue Probenschleife ein anderes Volumen hat als die Probenschleife, die Sie ausgebaut haben:
  - ♦ Prüfen Sie in der Konfiguration des Autosamplers auf der Seite Options, dass s das Volumen der Probenschleife korrekt angegeben ist (→ Seite 43). Ändern Sie die Einstellung, falls erforderlich. Übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Klicken Sie auf der Seite General auf die Schaltfläche Send configuration to module (→ Seite 45).
  - ♦ Prüfen Sie, dass eine geeignete Pufferschleife und Spritze installiert ist (→ Seite 17). Das Volumen der Probenschleife sollte mindestens dem Volumen der Spritze entsprechen.
- 12. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

# 7.7 Nadelsitz

#### 7.7.1 Wechseln des Nadelsitzes

Beschreibung	BestNr.
WPS-3000SL Analytical, WPS-3000TBSL Analytical, WPS-3000RS, WPS-3000TXRS: Nadelsitz (2 Stück) Servicekit Nadelsitz (1 Nadelsitz, 1 Halteschraube)	6820.0047A 6820.0038A
<i>WPS-3000TBRS:</i> Nadelsitz (2 Stück) Servicekit Nadelsitz (1 Nadelsitz, 1 Halteschraube)	6841.0047 6841.0038
WPS-3000SL Semiprep: Servicekit Nadelsitz (1 Nadelsitz, 1 Halteschraube)	6820.0043

- Bewegen Sie die Probennadel in die geeignete Position f
  ür den Wechsel des Nadelsitzes (Probennadel oben und Nadelarm ganz links).
   Rufen Sie am Ger
  ät das Men
  ü Maintenance auf und w
  ählen Sie Needle Position und Maintain— oder —geben Sie in Chromeleon das Kommando InitiateChangeNeedle.
- 2. Klappen Sie den Frontdeckel nach oben auf.
- 3. Entfernen Sie die Nadelsitzkapillare unten an der Waschport/Nadelsitz-Einheit mit einem Maulschlüssel (Größe 1/4", im Zubehör enthalten).



Abb. 62: Entfernen der Nadelsitzkapillare

4. Lösen Sie die Halteschraube mit einem Steckschlüssel (Größe 1/4").



Abb. 63: Entfernen der Halteschraube

- Entfernen Sie den Nadelsitz. 5.
- 6. Geben Sie einen Tropfen Flüssigkeit (Isopropanol oder Wasser) in die Halteschraube.



Abb. 64: Fixieren des Nadelsitzes

Setzen Sie den neuen Nadelsitz mit einer Pinzette in die Halteschraube ein. Achten Sie 7. auf die Ausrichtung: Der Konus muss in Richtung der kleinen Öffnung der Halteschraube zeigen ( $\rightarrow$  Abb. 65 und Abb. 66).



Beim Einsetzen in die Halteschraube muss der Konus des Nadelsitzes in Richtung der kleinen Öffnung der Schraube zeigen.

Abb. 65: Ausrichtung des Nadelsitzes



Setzen Sie den Nadelsitz so in die Halteschraube ein, dass der Konus in diese Richtung zeigt.

Abb. 66: Einsetzen des Nadelsitzes

8. Setzen Sie die Halteschraube auf und ziehen Sie die Schraube fest. Die Halteschraube muss bündig mit der Oberkante der Waschport/Nadelsitz-Einheit abschließen.



Abb. 67: Einsetzen des Nadelsitzes

Die Halteschraube muss mit dieser Kante bündig abschließen.

Nadelsitzkapillare

- 9. Schrauben Sie die Nadelsitzkapillare wieder an der Waschport/Nadelsitz-Einheit fest.
- 10. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder* geben Sie in Chromeleon das Kommando **SelfTest**.
- 11. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl NeedleSeatChanged.
- Empfehlung Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch (→ Seite 90).
- **I** Hinweis: Luftblasen in der Nadelsitzverbindung können die Verschleppung zeitweise erhöhen. Um die Luftblasen zu entfernen, schalten Sie den Fluss der Pumpe aus, schalten Sie das Injektionsventil in die Position **Inject** und lösen Sie die Nadelsitzkapillare an der Nadelsitz-Einheit mit einer 1/2-Umdrehung. Entlüften Sie das System über die Pumpe des UltiMate 3000-Systems. Damit der Leaksensor nicht anspricht, nehmen Sie die Flüssigkeit, die aus der Nadelsitzkapillare austritt, mit einem Tuch auf. Wenn der Entlüftungsvorgang abgeschlossen ist, stellen Sie den Pumpenfluss erneut aus und ziehen Sie die Nadelsitzkapillare wieder fest.

#### 7.7.2 Wechseln der Nadelsitzkapillare

Die Schritte zum Wechseln der Nadelsitzkapillare hängen vom Autosamplermodell ab. Folgen Sie den Schritten für Ihren Autosampler:

- WPS-3000SL Analytical, WPS-3000TBSL Analytical, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS: Folgen Sie den Schritten im Kapitel 7.7.2.1.
- WPS-3000SL Semiprep: Folgen Sie den Schritten im Kapitel 7.7.2.2 ( $\rightarrow$  Seite 144).

# 7.7.2.1 WPS-3000SL Analytical, WPS-3000TBSL Analytical, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS

Beschreibung	BestNr.
WPS-3000SL Analytical und WPS-3000RS: Nadelsitzkapillare, 3,1 µL, SST, Viper	6820.2464
WPS-3000TBSL Analytical: Nadelsitzkapillare, 3,1 µL, PEEK, Viper	6827.2408
WPS-3000TBRS: Nadelsitzkapillare, 1 μL, MP35N, Viper	6841.2472
WPS-3000TXRS: Nadelsitzkapillare, 1 μL, SST, Viper	6820.2472

- Bewegen Sie die Probennadel in die geeignete Position f
  ür den Wechsel des Nadelsitzkapillare (Probennadel oben und Nadelarm ganz links).
   Rufen Sie am Ger
  ät das Men
  ü Maintenance auf und w
  ählen Sie Needle Position und Maintain — oder — geben Sie in Chromeleon das Kommando InitiateChangeNeedle.
- 2. Klappen Sie den Frontdeckel nach oben auf.
- 3. Entfernen Sie die Nadelsitzkapillare unten an der Waschport/Nadelsitz-Einheit mit einem Maulschlüssel (Größe 1/4", im Zubehör enthalten).



Abb. 68: Entfernen der Nadelsitzkapillare

- 4. Entfernen Sie die Nadelsitzkapillare am Injektionsventil.
- 5. Führen Sie die neue Nadelsitzkapillare von unten in die Waschport/Nadelsitz-Einheit ein.



Abb. 69: Einführen der Nadelsitzkapillare

6. Ziehen Sie die Nadelsitzkapillare mit dem Maulschlüssel (Größe 1/4") fest.

Waschport/Nadelsitz-Einheit

Nadelsitzkapillare



- 7. Schließen Sie die neue Nadelsitzkapillare am Injektionsventil (Port 3) an.
- 8. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder*—geben Sie in Chromeleon das Kommando **SelfTest**.
- 9. Prüfen Sie die Verbindungen optisch auf Dichtigkeit.
- 10. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

Hinweis: Luftblasen in der Nadelsitzverbindung können die Verschleppung zeitweise erhöhen. Um die Luftblasen zu entfernen, schalten Sie den Fluss der Pumpe aus, schalten Sie das Injektionsventil in die Position Inject und lösen Sie die Nadelsitzkapillare an der Nadelsitz-Einheit mit einer 1/2-Umdrehung. Entlüften Sie das System über die Pumpe des UltiMate 3000-Systems. Damit der Leaksensor nicht anspricht, nehmen Sie die Flüssigkeit, die aus der Nadelsitzkapillare austritt, mit einem Tuch auf. Wenn der Entlüftungsvorgang abgeschlossen ist, stellen Sie den Pumpenfluss erneut aus und ziehen Sie die Nadelsitzkapillare wieder fest.

#### 7.7.2.2 WPS-3000SL Semiprep

Beschreibung	BestNr.
WPS-3000SL Semiprep: Nadelsitzkapillare, 24 μL, SST	6820.2409

1. Bewegen Sie die Probennadel in die geeignete Position für den Wechsel des Nadelsitzkapillare (Probennadel oben und Nadelarm ganz links).

Rufen Sie am Gerät das Menü **Maintenance** auf und wählen Sie **Needle Position** und **Maintain**— *oder* — geben Sie in Chromeleon das Kommando **InitiateChangeNeedle**.

- 2. Klappen Sie den Frontdeckel nach oben auf.
- 3. Entfernen Sie die Nadelsitzkapillare unten an der Waschport/Nadelsitz-Einheit mit einem Maulschlüssel (Größe 1/4", im Zubehör enthalten).

Entfernen Sie die Nadelsitzkapillare an der Nadelsitz-Einheit.



Abb. 71: Entfernen der Nadelsitzkapillare

- 4. Entfernen Sie die Nadelsitzkapillare am Injektionsventil.
- 5. Schieben Sie das Fitting und den Schneidring auf die neue Nadelsitzkapillare (Waschport/Nadelsitz-Ende). An den beiden Enden der Nadelsitzkapillare werden unterschiedliche Fittingverbindungen verwendet. Diese dürfen nicht vertauscht werden.



Abb. 72: Nadelsitzkapillare mit Fitting und Schneidring

6. Führen Sie die Nadelsitzkapillare von unten in die Waschport/Nadelsitz-Einheit ein.



Abb. 73: Führen der Nadelsitzkapillare

7. Schrauben Sie die Fittingschraube der Nadelsitzkapillare an der Waschport/Nadelsitz-Einheit zunächst lose an. (Die Nadelsitzkapillare kann dabei gegebenenfalls entnommen werden; sie wurde im vorherigen Schritt nur zum leichteren Einführen der Fittingverbindung benötigt.)



Abb. 74: Anziehen der Fittingschraube

8. Richten Sie die Nadelsitzkapillare aus (→ Abb. 75). Drücken Sie die Kapillare mit einer Pinzette fest in die Fittingverbindung.



Abb. 75: Ausrichten der Nadelsitzkapillare

 Ziehen Sie das Fitting der Nadelsitzkapillare mit dem Maulschlüssel fest (Größe 1/4"). Die Kapillare dreht sich gegebenenfalls am Schluss etwas mit. Zum Schutz der Kapillare sollten Sie dies nicht verhindern.



Abb. 76: Festziehen der Nadelsitzkapillare

10. Schieben Sie das Fitting und den Schneidring auf das andere Ende der Nadelsitzkapillare und führen Sie die Kapillare vorsichtig zum Injektionsventil.



Schieben Sie dieses Fitting auf das Injektionsventil-Ende der Kapillare.



Abb. 77: Nadelsitzkapillare mit Fittingverbindung

11. Schließen Sie die Nadelsitzkapillare am Injektionsventil (Port 3) an. Drücken Sie die Kapillare mit einer Pinzette bis zum Anschlag in den Port. Ziehen Sie die Kapillare mit dem Maulschlüssel fest (Größe 1/4").



Abb. 78: Befestigen der Nadelsitzkapillare am Injektionsventil

- 12. Fahren Sie den Nadelarm und die Probennadel in die korrekte Ausgangsposition zurück. Rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Self test**—*oder* geben Sie in Chromeleon das Kommando **SelfTest**.
- 13. Prüfen Sie die Verbindungen optisch auf Dichtigkeit.
- 14. Empfehlung

Testen Sie den Autosampler auf Dichtigkeit. Führen Sie in Chromeleon den Leak Test durch ( $\rightarrow$  Seite 90).

**I** Hinweis: Luftblasen in der Nadelsitzverbindung können die Verschleppung zeitweise erhöhen. Um die Luftblasen zu entfernen, schalten Sie den Fluss der Pumpe aus, schalten Sie das Injektionsventil in die Position **Inject** und lösen Sie die Nadelsitzkapillare an der Nadelsitz-Einheit mit einer 1/2-Umdrehung. Entlüften Sie das System über die Pumpe des UltiMate 3000-Systems. Damit der Leaksensor nicht anspricht, nehmen Sie die Flüssigkeit, die aus der Nadelsitzkapillare austritt, mit einem Tuch auf. Wenn der Entlüftungsvorgang abgeschlossen ist, stellen Sie den Pumpenfluss erneut aus und ziehen Sie die Nadelsitzkapillare wieder fest.

## 7.8 Injektionsventil

Der Autosampler ist mit einem Injektionsventil (2 Positionen, 6 Ports) ausgestattet ( $\rightarrow$  Seite 27). Beachten Sie die folgenden Hinweise für einen optimalen Betrieb des Injektionsventils:

- Das Ventil sollte nicht mit Lösungen verwendet werden, die konzentrierte Ameisensäure (Methansäure) enthalten.
- Um eine Beeinträchtigung der Dichtflächen des Ventils zu vermeiden, sollten Sie das Ventil nicht bei Temperaturen unter 0 °C mit Wasser in der Fluidik betreiben, lagern oder versenden.
- Verwenden Sie für die Anschlüsse am Injektionsventil möglichst Viper-Kapillarverbindungen bzw. ausschließlich die Schneidringe und Fittingschrauben aus dem Gerätezubehör oder Original-Dionex-Ersatzteile.



Abb. 79: Viper-Fttingverbindungen und andere Ventilschneidringe (Beispiele)

Der Wartungsaufwand für das Ventil ist sehr gering. In den meisten Fällen, reicht es aus, das Ventil gründlich mit einem geeigneten Lösungsmittel zu spülen. Dabei ist die Wahl des Mittels abhängig von der Art der Probe und der verwendeten mobilen Phase.

Die Dichtungen (Rotor Seal und Stator Face Seal) sind Verschleißteile und sollten nach Bedarf gewechselt werden ( $\rightarrow$  Seite 155).

#### 7.8.1 Prüfen des Injektionsventils auf Undichtigkeit

Auf dem Statorring des Injektionsventils befinden sich mehrere Auslassöffnungen. Tritt an einer oder mehreren Auslassöffnungen Flüssigkeit aus, kann dies auf eine mögliche Undichtigkeit des Ventils hindeuten.



Abb. 80: Auslassöffnungen am Injektionsventil (Beispiel)

Folgende Abhilfemaßnahmen sind möglich:

- Ersetzen Sie das Rotor Seal bzw. das Rotor Seal und Stator Seal ( $\rightarrow$  Seite 155).
- Ersetzen Sie den Stator ( $\rightarrow$  Seite 154).
- Ersetzen Sie den Ventil-Pod ( $\rightarrow$  Seite 150).

#### 7.8.2 Wechseln des Ventil-Pods

Beschreibung	BestNr.
<ul> <li>WPS-3000SL Analytical und WPS-3000SL Semiprep:</li> <li>Ventil-Pod (Injektionsventil)</li> <li>Der Pod ist geeignet für einen Druck &lt; 62 MPa (9000 psi)</li> <li>Das Rotor-Seal im Ventil-Pod ist geeignet für pH-Werte 0 - 14.</li> </ul>	6822.0051
<i>WPS-3000TBSL Analytical:</i> Ventil-Pod (Injektionsventil) Der Pod ist geeignet für einen Druck < 34 MPa (5000 psi) Das Rotor-Seal im Ventil-Pod ist geeignet für pH-Werte 0 - 14.	6723.9013
<i>WPS-3000RS:</i> Ventil-Pod (Injektionsventil) Der Pod ist geeignet für einen Druck <103 MPa (15000 psi) Das Rotor-Seal im Ventil-Pod ist für pH-Werte 0 - 14 geeignet.	6840.0013
<i>WPS-3000TBRS:</i> Ventil-Pod (Injektionsventil) Der Pod ist geeignet für einen Druck <103 MPa (15000 psi) Das Rotor-Seal im Ventil-Pod ist für pH-Werte 0 - 14 geeignet.	6841.0013
<i>WPS-3000TXRS:</i> Ventil-Pod (Injektionsventil) Der Pod ist geeignet für einen Druck <125 MPa (18130 psi) Das Rotor-Seal im Ventil-Pod ist für pH-Werte 0-10 geeignet.	6843.0013

 Schalten Sie das Injektionsventil in die Position Inject.
 Rufen Sie am Gerät das Menü Maintenance auf und wählen Sie Inject valve und Inject— oder — geben Sie in Chromeleon das Kommando InjectValveToInject.

- 2. Lösen Sie alle Verbindungen am Injektionsventil.
- 3. Folgen Sie der Beschreibung für Ihren Autosampler:
  - WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000TBRS, WPS-3000RS und WPS-3000TXRS: Folgen Sie der Beschreibung auf Seite 151.
  - WPS-3000TBSL Analytical: Folgen Sie der Beschreibung auf Seite 152.

# WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS

Dieser Abschnitt gilt nicht für den WPS-3000TBSL Analytical. Für diesen Autosampler, siehe ab Seite 152.

4. Lösen Sie die Überwurfmutter ("Spanner"). Lösen Sie die Überwurfmutter nur mit der Hand (verwenden Sie kein Werkzeug).



Abb. 81: Lösen der Überwurfmutter

- 5. Ziehen Sie den Pod nach unten ab.
- 6. Richten Sie den Positionierstift im neuen Pod mit der entsprechenden Aussparung im Ventilkörper aus. Schieben Sie den Pod vollständig in den Ventilkörper hinein.



Abb. 82: Ausrichten des Positionierstifts

Wenn dies nicht möglich ist, folgen Sie den Schritten unter Troubleshooting-Tip (siehe unten).

- 7. Ziehen Sie die Überwurfmutter manuell (ohne Werkzeug) handfest an. Wird die Überwurfmutter zu stark anzogen, lässt sich das Ventil eventuell nicht mehr schalten. Beim Versuch, das Ventil zu schalten, erscheint eine Meldung, z.B. Motor switching valve malfunction.
- 8. Mit einem Wechsel des Ventil-Pods wird automatisch auch das Rotor Seal gewechselt. Aktualisieren Sie daher in Chromeleon die Serviceinformationen für das Rotor Seal. Geben Sie den Befehl **RotorSealChanged**.

#### **Troubleshooting-Tip**

Wenn Sie den Pod nicht vollständig in den Ventilkörper schieben können, stimmt die Stellung der Anschlusskupplung nicht mit der aktuellen Position des Antriebs überein. Gehen Sie folgendermaßen vor:

- 1. Schalten Sie das Ventil in die Position Load und versuchen Sie es erneut.
- 2. Wenn Sie den Pod auch jetzt nicht vollständig in den Ventilkörper schieben können, drehen Sie den Pod zunächst im Ventilkörper, bis die Anschlusskupplung im Antrieb einrastet. Richten Sie dann durch Weiterdrehen den Positionierstift mit der Aussparung aus und schieben Sie den Pod vollständig in den Ventilkörper hinein.

#### WPS-3000TBSL Analytical

4. Lösen Sie die Überwurfmutter ("Spanner"). Lösen Sie die Überwurfmutter nur mit der Hand (verwenden Sie kein Werkzeug).



Abb. 83: Lösen der Überwurfmutter

- 5. Ziehen Sie den Pod nach unten ab.
- 6. Schieben Sie den neuen Pod vollständig in den Ventilkörper hinein.
   Achten Sie auf die korrekte Ausrichtung. Richten Sie die *breite* Vertiefung im Pod
   (→ Abb. 84) mit der entsprechenden Aufnahme im Ventilkörper aus.



Aufnahme im Ventilkörper

Breite Vertiefung im Pod

Abb. 84: Einbau des Pods

Wenn Sie den Pod nicht richtig in den Ventilkörper schieben können, folgen Sie den Schritten unter Troubleshooting-Tip ( $\rightarrow$  Seite 153).

- 7. Ziehen Sie die Überwurfmutter manuell (ohne Werkzeug) handfest an. Wird die Überwurfmutter zu stark anzogen, lässt sich das Ventil eventuell nicht mehr schalten. Beim Versuch, das Ventil zu schalten, erscheint eine Meldung, z.B. Motor switching valve malfunction.
- 8. Mit einem Wechsel des Ventil-Pods wird automatisch auch das Rotor Seal gewechselt. Aktualisieren Sie daher in Chromeleon die Serviceinformationen für das Rotor Seal. Geben Sie den Befehl **RotorSealChanged**.

#### **Troubleshooting-Tip**

Wenn Sie den Pod nicht richtig in den Ventilkörper schieben können, stimmt die Stellung der Anschlusskupplung nicht mit der aktuellen Position des Antriebs überein. Gehen Sie folgendermaßen vor:

- 1. Schalten Sie das Ventil in die Position Load und versuchen Sie es erneut.
- 2. Wenn Sie den Pod auch jetzt nicht vollständig in den Ventilkörper schieben können, ist die Anschlusskupplung verdreht. Drehen Sie die Anschlusskupplung manuell, bis Sie den Pod in den Ventilkörper schieben können.

#### 7.8.3 Wechseln des Ventilstators

Beschreibung	BestNr.
WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000RS: Stator (Injektionsventil)	6840.0011
WPS-3000TBSL Analytical: Stator (Injektionsventil)	6723.9011
WPS-3000TBRS: Stator (Injektionsventil)	6841.0011
WPS-3000TXRS:Stator (Injektionsventil)Empfehlung: Nach einem Tausch des Stators wird unter Umständen nicht immer dievolle Lebensdauer des Ventils erreicht. Thermo Fisher Scientific empfiehlt daher, denVentil-Pod zu tauschen ( $\rightarrow$ Seite 154).	6840.0011

- 1. Entfernen Sie den Ventil-Pod:
  - WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000TBRS, WPS-3000RS und WPS-3000TXRS: Folgen Sie der Beschreibung auf Seite 151.
  - WPS-3000TBSL Analytical: Folgen Sie der Beschreibung auf Seite 152.
- Entfernen Sie die drei Statorschrauben mit einem Innensechskantschlüssel (Größe 3/32"). Um ein Verkanten des Stators zu vermeiden, entfernen Sie die drei Schrauben gleichzeitig.



Statorschrauben

*Abb.* 85: *Statorschrauben (hier am Edelstahlventil)* 

- 3. Entfernen Sie den Stator vom Ventilkörper.
- 4. Setzen Sie den neuen Stator so auf den Ventilkörper auf, dass die Befestigungspins im Ventilkörper in die entsprechenden Bohrungen im Stator passen.
- 5. Ziehen Sie die Schrauben zunächst abwechselnd handfest an. Ziehen Sie die Schrauben dann fest, dabei sollte das Drehmoment 0,7 Nm nicht übersteigen.
- 6. Bauen Sie den Pod wieder ein.

#### 7.8.4 Rotor Seal und Stator Face Seal

Beschreibung	BestNr.
WPS-3000SL Analytical und WPS-3000SL Semiprep , WPS-3000RS und WPS-3000TBRS:	
Rotor Seal (geeignet für pH-Werte 0 - 14)	6840.0012
<i>WPS-3000TBSLAnalytical:</i> Rotor Seal und Stator Face Seal (für pH-Werte 0 – 14)	6722.9014
WPS-3000TXRS: Rotor Seal (geeignet für pH-Werte 0-10)	6843.0012
<i>Empfehlung</i> : Nach einem Tausch des Rotors Seals wird unter Umständen nicht immer die volle Lebensdauer des Ventils erreicht. Thermo Fisher Scientific empfiehlt daher, den Ventil-Pod zu tauschen ( $\rightarrow$ Seite 154).	

Folgen Sie den Schritten für Ihren Autosampler:

- WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS: Folgen Sie den Schritten in Kapitel 7.8.4.1 (→ Seite 156).
- WPS-3000TBSL Analytical: Folgen Sie den Schritten in Kapitel 7.8.4.2 (→ Seite 157).

Beachten Sie beim Wechseln der Dichtungen Folgendes:

- Um eine Beschädigung der Dichtflächen des Ventils zu vermeiden, verwenden Sie beim Entfernen und Einbauen einer Dichtung keine spitzen oder scharfkantigen Werkzeuge. Selbst kleinste Kratzer können die Dichtigkeit des Ventils beeinträchtigen.
- Achten Sie beim Einbau der neuen Dichtung darauf, dass keine Verunreinigungen an den Ventilteilen haften. Selbst kleinste Partikel können zu Schäden am Ventil führen und die Dichtigkeit beeinträchtigen.

# 7.8.4.1 WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS

Dieses Kapitel gilt nicht für den WPS-3000TBSL Analytical. Für diesen Autosampler, siehe Kapitel 7.8.4.2 ( $\rightarrow$  Seite 157).

- 1. Entfernen Sie den Ventil-Pod ( $\rightarrow$  Seite 151).
- 2. Entfernen Sie den Stator ( $\rightarrow$  Seite 154).
- Entfernen Sie das Rotor Seal. Beachten Sie dabei die Hinweise auf Seite 155.



Abb. 86: Ventilkörper ohne Rotor Seal und Stator

4. Setzen Sie das neue Rotor Seal so auf den Ventilkörper auf, dass die Gravuren im Seal zum Stator zeigen. Die drei Stifte im Ventilkörper passen nur in einer Richtung in die entsprechenden Löcher im Rotor Seal. Beachten Sie auch die Hinweise auf Seite 155.



Abb. 87: Rotor Seal

- 5. Bauen Sie den Stator wieder ein ( $\rightarrow$  Seite 154).
- 6. Bauen Sie den Pod wieder ein ( $\rightarrow$  Seite 150).
- 7. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl **RotorSealChanged**.

#### 7.8.4.2 WPS-3000TBSL Analytical

- 1. Entfernen Sie den Ventil-Pod ( $\rightarrow$  Seite 152).
- 2. Entfernen Sie den Stator ( $\rightarrow$  Seite 154).
- Entfernen Sie das Rotor Seal. Beachten Sie dabei die Hinweise auf Seite 155.
- 4. Setzen Sie das neue Rotor Seal so auf die Aufnahme, dass die Gravuren zum Stator zeigen. Die drei Stifte in der Aufnahme passen nur in einer Richtung in die entsprechenden Löcher im Rotor Seal (siehe Montageanleitung aus dem Kit). Beachten Sie auch die Hinweise auf Seite 155.
- 5. Entfernen Sie das alte Stator Face Seal aus dem Stator. Beachten Sie dabei die Hinweise auf Seite 155.
- 6. Setzen Sie das neue Stator Face Seal in den Stator ein. Die drei Stifte in der Dichtung passen nur in einer Richtung in die entsprechenden Löcher im Stator. Beachten Sie auch die Hinweise auf Seite 155.
- 7. Bauen Sie den Stator wieder ein ( $\rightarrow$  Seite 154).
- 8. Bauen Sie den Pod wieder ein ( $\rightarrow$  Seite 152).
- 9. Aktualisieren Sie die Serviceinformationen in Chromeleon. Geben Sie den Befehl **RotorSealChanged**.

# 7.9 Wechseln der Sicherungen

**Warnung:** Schalten Sie den Autosampler aus. Ziehen Sie den Netzstecker.

1. Hebeln Sie mit einem kleinen Schraubendreher den Sicherungshalteschlitten aus der Netzbuchse.



Abb. 88: Sicherungshalteschlitten

2. Wechseln Sie die Sicherungen.



**ng:** Setzen Sie immer zwei neue Sicherungen ein. Verwenden Sie nur die unten angegebenen Sicherungen.

Beschreibung	BestNr.
Autosampler ohne Probenthermostatisierung:	Enthalten in: WPS-Sicherungskit,
Sicherung, 2 A, träge, 5 x 20mm	BestNr. 6820.0026
Autosampler mit Probenthermostatisierung:	Informationen zum Inhalt des Kits finden
Sicherung, 4 A, träge, 5 x 20mm	Sie im Kapitel 10.3 (→ Seite 183).

- 3. Setzen Sie den Sicherungshalteschlitten wieder auf.
- 4. Stecken Sie das Netzkabel wieder an. Schalten Sie den Autosampler ein.

### 7.10 Aktualisieren der Autosampler-Firmware

Die aktuelle Firmware-Version ist bei Auslieferung des Autosamplers installiert. Die Autosampler-Firmware ist auch in Chromeleon enthalten.

Welche Firmware-Version im Autosampler installiert ist und welche in Chromeleon enthalten ist, stellen Sie so fest:

- Firmware-Version des Autosamplers
  - Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite ein. Auf dem Autosamplerdisplay erscheinen allgemeine Informationen zum Autosampler, einschließlich der Firmware-Version.
  - Rufen Sie am Autosamplerdisplay das Menü **Diagnostics** auf (→ Seite 79) und wählen Sie **Firmware version**.
- *Firmware-Version in Chromeleon* Navigieren Sie im Windows-Explorer zu der Datei **IQReport.log** im Ordner **IQ** Ihrer Chromeleon-Installation. Suchen Sie in der Datei nach UM3\_WPS.hex.

**i** Hinweis: Die Informationen zu den Firmware-Versionen erhalten Sie auch, wenn Sie die Autosampler-Firmware über Chromeleon aktualisieren (siehe unten).

Wenn eine neue Firmware-Version für den Autosampler verfügbar ist, wird diese zusammen mit der nächsten Service Release zu Chromeleon ausgeliefert und in den Release Notes beschrieben.

Wenn Sie die Service Release installieren, wird die neue Firmware *nicht* automatisch auf den Autosampler übertragen. Übertragen Sie die neue Firmware wie folgt:

Vorsicht: Damit die Aktualisierung erfolgreich durchgeführt werden kann, darf die Kommunikation zwischen Chromeleon und dem Autosampler während der Übertragung *keinesfalls* unterbrochen oder der Autosampler ausgeschaltet werden.

- 1. Vergewissern Sie sich, dass folgende Voraussetzungen erfüllt sind:
  - Der Autosampler ist in Chromeleon verbunden ("connected").
  - Der Chromeleon-Server befindet sich im Modus *running idle*. Es laufen keine Prozesse auf dem Chromeleon Server-PC oder in Chromeleon.
- 2. Starten Sie das Programm Server Configuration ( $\rightarrow$  Seite 39).
- 3. Markieren Sie den Autosampler in der Zeitbasis mit einem Rechtsklick und wählen Sie im Menü den Punkt **Properties**.

- Klicken Sie auf der Registerkarte General (→ Seite 40) auf die Schaltfläche Firmware Download. Eine Meldung informiert Sie über die aktuell im Autosampler installierte Firmware-Version und gibt an, welche Version bei einem Download von Chromeleon auf den Autosampler übertragen wird.
  - **i** Hinweis: Wenn die Firmware im Autosampler eine neuere Version ist als in Chromeleon, sollten Sie die Firmware-Version aus Chromeleon *nicht* auf den Autosampler übertragen. Ältere Firmware-Versionen sind gegebenenfalls nicht mit neuerer Hardware kompatibel.
- 5. Klicken Sie **Yes**, wenn Sie die Übertragung starten möchten. (Klicken Sie **No**, wenn Sie die Aktualisierung nicht durchführen möchten.)

Die Übertragung kann einige Minuten dauern. Sie ist abgeschlossen, wenn in der Serverkonfiguration im Fenster **Messages Server** die Meldung **Download completed successfully** erscheint. Die Meldung erscheint auch im Chromeleon Audit Trail.

Wird der Download nicht erfolgreich abgeschlossen, erscheinen entsprechende Meldungen im Audit Trail. Schalten Sie den Autosampler in diesem Fall aus und wieder ein. Führen Sie den Download wie oben beschrieben erneut durch. Führt dies nicht zum Erfolg, wenden Sie sich an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.

**i** Hinweis: Nach Abschluss des Downloads von Chromeleon wird gegebenenfalls noch ein interner Download am Gerät durchgeführt. Ist dies der Fall, können Sie den Fortschritt am Gerätedisplay verfolgen. Schalten Sie den Autosampler keinesfalls aus, bis der interne Download abgeschlossen ist.

# 8 Optimieren für spezielle Anwendungen

Um den WPS-3000RS für Anwendungen in der Mikro-HPLC zu optimieren, benötigen Sie:

Beschreibung	BestNr.
Probenschleife Mikro, 25 $\mu$ L (V = 40 $\mu$ L), SST, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen.	6820.2452

- 1. Wechseln Sie die Probenschleife ( $\rightarrow$  Seite 131).
- 2. Ändern Sie die Konfiguration des Autosamplers in Chromeleon:
  - a) Öffnen Sie das Dialogfenster **Properties** für den Autosampler ( $\rightarrow$  Seite 45).
  - b) Klicken Sie auf der Seite **Options** das Kontrollkästchen **Micro** an ( $\rightarrow$  Seite 43).
- 3. Übertragen Sie die geänderte Konfiguration von Chromeleon zum Autosampler. Klicken Sie **Send configuration to module** auf der Seite **General** (→ Seite 45).

# 9 Technische Daten

Alle technische Daten: Stand September 2013 - Änderungen vorbehalten

# 9.1 WPS-3000SL Analytical und WPS-3000SL Semiprep

	WPS-3000(T)SL Analytical	WPS-3000(T)SL Semiprep
Probenkapazität	Abhängig von der Drehtellerkonfiguration:	
Fläschchen	<ul> <li>216 x 0,3-mL-Fläschchen, 120 x 1,1-mL konische Fläschchen,</li> <li>216 x 1,2-mL-Fläschchen, 120 x 1,8- mL/2,0-mL-Fläschchen,</li> <li>66 x 4-mL-Fläschchen, 30 x 10- mL-Fläschchen bzw.</li> <li>120 x 0,5-mL oder 1,5-mL Eppendorf Fläschchen (offen, ohne Kappe)</li> <li>+ 15 x 10 mL Probenfläschchen</li> </ul>	
Well Plates	3 x 96 bzw. 3 x 384 Mikrotiterplatte 3 x 24 Deep-Well-Plates, 3 x 384 Lo + 15 x 10 mL Probenfläschchen	en oder Deep-Well-Plates, ow-Well-PCR-Plates
Injektionsmethode	Inline-Split-Loop	
<b>Injektionsvolumen</b> Einstellbar (empfohlen) Optional	0,01 – 100 μL (1 – 100 μL) 0,2 – 25 μL, 1,5 – 250 μL, 1,5 – 500 μL	0,01 – 2500 μL (100 – 2000 μL) 10 – 1000 μL
Genauigkeit Injektionsvolumen	Typisch $\pm 0,5\%$ bei 50 und 90 $\mu$ L	Typisch ±1% bei 2000 µL
Erforderliche Probenmenge (Minimum)	1 μL von 5 μL (250-μL-Fläschchen, konisch)	Entfällt
Präzision Injektionsvolumen	<0,25% RSD bei 5 µL (typisch <0,15% RSD), Koffein in Wasser	<0,3% RSD bei 100 µL (typisch <0,15% RSD), Koffein in Wasser
Linearität	Korr. koeff. >0,9999, RSD <0.5% bei 5-90 μL, Koffein in Wasser	Korr. koeff. >0,9995, RSD <1% bei 100-2000 μL, Koffein in Wasser
Waschen der Nadel	Aktives Waschen der Probennadel von außen	
Verschleppung	<0,004% für Koffein mit externem Waschen bei 7.5 MPa	<0,004% für Koffein mit externem Waschen bei 7.5 MPa
Injektionszykluszeit	<15 s für 5 µL	<20 s für 100 μL
<b>WPS-3000TSL:</b> Probentemperierung Genauigkeit Probentemperatur	4 bis 45 °C oder 22 °C unterhalb der Umgebungstemperatur ±2 °C bei einem Sollwert von 10 °C bei Umgebungstemperaturen von <25 °C und <50 % relativer Luftfeuchtigkeit	
Gewicht	19 kg (WPS-3000SL) 24 kg (WPS-3000TSL)	

	WPS-3000(T)SL Analytical	WPS-3000(T)SL Semiprep
Abmessungen (H x B x T)	$36 \times 42 \times 51 \text{ cm}$	
PC-Anschluss	Alle Funktionen und Parameter software-gesteuert über USB 1.1 Integrierter USB-Hub mit drei USB 1.1-Schnittstellen	
E/A-Schnittstellen	4 digitale Eingänge, 4 programmiert	oare Relaisausgänge
Sicherheitsmerkmale	Leaksensor, automatische Platten-/P interne Überwachung aller mechanis WPS-3000TSL: Überwachung der F Temperatursensoren	robenfläschchenerkennung, schen Operationen funktion und Kühlleistung über 3
Eingabe/Anzeige	LCD zur Anzeige von Systemparametern, Standby-Taste 3 LEDs (Power, Connected und Status) zur Statusangabe 4 Funktionstasten zum Betrieb während der Erstinstallation und zur Wartung	
GLP	Unter Chromeleon: Automatische G Routinen) und Überwachung der Sy Alle Systemparameter werden im Cl	erätequalifikation (AutoQ <sup>™</sup> - stem Wellness. nromeleon Audit Trail protokolliert.
Medienberührte Teile		
Flusspfad Probe	PEEK <sup>™</sup> , Edelstahl	
Flusspfad Waschflüssigkeit	PEEK, Edelstahl, PCTFE, Fused Sil	ica, ETFE, ECTFE, PTFE, Glas
Leistungsaufnahme	100–120 V, 60 Hz; 200–240 V, 50 F max. 150 VA (WPS-3000SL) bzw. 1	Hz; nax. 320 VA (WPS-3000TSL)
Emissionsschalldruckpegel	Typisch $< 65 \text{ dB}(A)$ in 1 m Abstance	1
Umgebungsbedingungen	Verwendungsbereich: Innenraum Temperaturbereich: 10 °C bis 35 Luftfeuchtigkeit: 80% relative Überspannungskat.: II Verschmutzungsgrad: 2	°C Feuchte, nicht kondensierend

	WPS-3000TBSL Analytical
Probenkapazität	Abhängig von der Drehtellerkonfiguration:
Fläschchen	<ul> <li>216 x 0,3-mL-Fläschchen, 120 x 1,1-mL konische Fläschchen,</li> <li>216 x 1,2-mL-Fläschchen, 120 x 1,8- mL/2,0-mL-Fläschchen,</li> <li>66 x 4-mL-Fläschchen, 30 x 10- mL-Fläschchen bzw.</li> <li>120 x 0,5-mL oder 1,5-mL Eppendorf Fläschchen (offen, ohne Kappe)</li> <li>+ 15 x 10 mL Probenfläschchen</li> </ul>
Well Plates	3 x 96 bzw. 3 x 384 Mikrotiterplatten oder Deep-Well-Plates, 3 x 24 Deep-Well-Plates, 3 x 384 Low-Well-PCR-Plates + 15 x 10 mL Probenfläschchen
Injektionsmethode	Inline-Split-Loop
Injektionsvolumen Einstellbar (empfohlen)	0,01 – 100 μL (1 – 100 μL)
Genauigkeit Injektionsvolumen	Typisch $\pm 0.5\%$ bei 50 und 90 $\mu$ L
Erforderliche Probenmenge (Minimum)	1 μL von 5 μL (250-μL-Fläschchen, konisch)
Präzision Injektionsvolumen	<0,25% RSD bei 5 µL (typisch <0,15% RSD), Koffein in Wasser
Linearität	Korr. koeff. >0,9999, RSD <0.5% bei 5-90 μL, Koffein in Wasser
Waschen der Nadel	Aktives Waschen der Probennadel von außen
Verschleppung	<0,004% für Koffein mit externem Waschen bei 7.5 MPa
Injektionszykluszeit	<15 s für 5 μL
Probentemperierung	4 bis 45 °C oder 22 °C unterhalb der Umgebungstemperatur
Genauigkeit Probentemperatur	$\pm 2$ °C bei einem Sollwert von 10 °C bei Umgebungstemperaturen von $\leq 25$ °C und $\leq 50$ % relativer Luftfeuchtigkeit
Gewicht	24 kg
Abmessungen (H x B x T)	$36 \times 42 \times 51$ cm
PC-Anschluss	Alle Funktionen und Parameter software-gesteuert über USB 1.1 Integrierter USB-Hub mit drei USB 1.1-Schnittstellen
E/A-Schnittstellen	4 digitale Eingänge, 4 programmierbare Relaisausgänge
Sicherheitsmerkmale	Leaksensor, automatische Platten-/Probenfläschchenerkennung, interne Überwachung aller mechanischen Operationen Überwachung der Funktion und Kühlleistung über 3 Temperatursensoren
Eingabe/Anzeige	LCD zur Anzeige von Systemparametern, Standby-Taste 3 LEDs (Power, Connected und Status) zur Statusangabe 4 Funktionstasten zum Betrieb während der Erstinstallation und zur Wartung

	WPS-3000TBSL Analytical
GLP	Unter Chromeleon: Automatische Gerätequalifikation (AutoQ <sup>™</sup> - Routinen) und Überwachung der System Wellness. Alle Systemparameter werden im Chromeleon Audit Trail protokolliert.
Medienberührte Teile	
Flusspfad Probe	PEEK™, Edelstahl
Flusspfad Waschflüssigkeit	PEEK, Edelstahl, PCTFE, Fused Silica, ETFE, ECTFE, PTFE, Glas
Leistungsaufnahme	100–120 V, 60 Hz; 200–240 V, 50 Hz; max. 320 VA
Emissionsschalldruckpegel	Typisch $< 65 \text{ dB}(A)$ in 1 m Abstand
Umgebungsbedingungen	Verwendungsbereich:InnenraumTemperaturbereich:10 °C bis 35 °CLuftfeuchtigkeit:80% relative Feuchte, nicht kondensierendÜberspannungskat.:IIVerschmutzungsgrad:2

## 9.3 WPS-3000RS

	WPS-3000(T)RS
Probenkapazität	Abhängig von der Drehtellerkonfiguration:
Fläschchen	<ul> <li>216 x 0,3-mL-Fläschchen, 120 x 1,1-mL konische Fläschchen,</li> <li>216 x 1,2-mL-Fläschchen, 120 x 1,8- mL/2,0-mL-Fläschchen,</li> <li>66 x 4-mL-Fläschchen, 30 x 10- mL-Fläschchen bzw.</li> <li>120 x 0,5-mL oder 1,5-mL Eppendorf Fläschchen (offen, ohne Kappe)</li> <li>+ 15 x 10 mL Probenfläschchen</li> </ul>
Well Plates	3 x 96 bzw. 3 x 384 Mikrotiterplatten oder Deep-Well-Plates, 3 x 24 Deep-Well-Plates, 3 x 384 Low-Well-PCR-Plates + 15 x 10 mL Probenfläschchen
Injektionsmethode	Inline-Split-Loop
<b>Injektionsvolumen</b> Einstellbar (empfohlen) Optional	0,01 – 100 μL (1 – 100 μL) 0,2 – 25 μL, 1,5 – 250 μL, 1,5 – 500 μL
Genauigkeit Injektionsvolumen	Typisch $\pm 0.5\%$ bei 20 µL
Erforderliche Probenmenge (Minimum)	1 μL von 5 μL (250-μL-Fläschchen, konisch)
Präzision Injektionsvolumen	${<}0,\!25\%$ RSD bei 5 $\mu L$ (typisch ${<}0,\!15\%$ RSD), Koffein in Wasser
Linearität	Korr. koeff. >0,9999, RSD <0.5% bei 5-90 $\mu$ L, Koffein in Wasser
Waschen der Nadel	Aktives Waschen der Probennadel von außen
Verschleppung	<0,004% für Koffein mit externem Waschen bei 20 MPa
Injektionszykluszeit	<15 s für 5 μL
WPS-3000TRS	
Probentemperierung	4 bis 45 °C oder 22 °C unterhalb der Umgebungstemperatur
Genauigkeit Probentemperatur	$\pm 2$ °C bei einem Sollwert von 10 °C bei Umgebungstemperaturen von $\leq 25$ °C und $\leq 50$ % relativer Luftfeuchtigkeit
Gewicht	19 kg (WPS-3000RS) 24 kg (WPS-300TRS)
Abmessungen (H x B x T)	$36 \times 42 \times 51$ cm
PC-Anschluss	Alle Funktionen und Parameter software-gesteuert über USB 1.1 Integrierter USB-Hub mit drei USB 1.1-Schnittstellen
E/A-Schnittstellen	4 digitale Eingänge, 4 programmierbare Relaisausgänge
Sicherheitsmerkmale	Leaksensor, automatische Platten-/Probenfläschchenerkennung, interne Überwachung aller mechanischen Operationen WPS-3000TRS: Überwachung der Funktion und Kühlleistung über 3 Temperatursensoren
Eingabe/Anzeige	LCD zur Anzeige von Systemparametern, Standby-Taste 3 LEDs (Power, Connected und Status) zur Statusangabe 4 Funktionstasten zum Betrieb während der Erstinstallation und zur Wartung

	WPS-3000(T)RS	
GLP	Unter Chromeleon: Au Routinen) und Überwa Alle Systemparameter protokolliert.	tomatische Gerätequalifikation (AutoQ <sup>™</sup> - chung der System Wellness. werden im Chromeleon Audit Trail
Medienberührte Teile		
Flusspfad Probe	PEEK™, Edelstahl	
Flusspfad Waschflüssigkeit	PEEK, Edelstahl, PCT	FE, Fused Silica, ETFE, ECTFE, PTFE, Glas
Leistungsaufnahme	100–120 V, 60 Hz; 200 max. 150 VA (WPS-30	0–240 V, 50 Hz; 000RS) bzw. max. 320 VA (WPS-3000TRS)
Emissionsschalldruckpegel	Typisch $< 65 \text{ dB}(A)$ in	n 1 m Abstand
Umgebungsbedingungen	Verwendungsbereich: Temperaturbereich: Luftfeuchtigkeit: Überspannungskat.: Verschmutzungsgrad:	Innenraum 10 °C bis 35 °C 80% relative Feuchte, nicht kondensierend II 2

## 9.4 WPS-3000TBRS

	WPS-3000TBRS
Probenkapazität	Abhängig von der Drehtellerkonfiguration:
Fläschchen	<ul> <li>216 x 0,3-mL-Fläschchen, 120 x 1,1-mL konische Fläschchen,</li> <li>216 x 1,2-mL-Fläschchen, 120 x 1,8- mL/2,0-mL-Fläschchen,</li> <li>66 x 4-mL-Fläschchen, 30 x 10- mL-Fläschchen bzw.</li> <li>120 x 0,5-mL oder 1,5-mL Eppendorf Fläschchen (offen, ohne Kappe)</li> <li>+ 15 x 10 mL Probenfläschchen</li> </ul>
Well Plates	3 x 96 bzw. 3 x 384 Mikrotiterplatten oder Deep-Well-Plates, 3 x 24 Deep-Well-Plates, 3 x 384 Low-Well-PCR-Plates + 15 x 10 mL Probenfläschehen
Injektionsmethode	Inline-Split-Loop
Injektionsvolumen Einstellbar (empfohlen) Optional	0,01 - 25 μL (0,2 – 25 μL) 1 – 100 μL, 1,5 – 250 μL, 1,5 – 500 μL
Genauigkeit Injektionsvolumen	Typisch ±0,5% bei 20 μL
Erforderliche Probenmenge (Minimum)	1 μL von 5 μL (250-μL-Fläschchen, konisch)
Präzision Injektionsvolumen	<0,3% RSD bei 2 $\mu$ L (typisch <0,15% RSD), Koffein in Wasser
Linearität	Korr. koeff. >0,9999, RSD <0.5% bei 1-20 $\mu$ L, Koffein in Wasser
Waschen der Nadel	Aktives Waschen der Probennadel von außen
Verschleppung	<0,004% für Koffein mit externem Waschen bei 20 MPa
Injektionszykluszeit	<20 s für 5 μL
Probentemperierung	4 bis 45 °C oder 22 °C unterhalb der Umgebungstemperatur
Genauigkeit Probentemperatur	$\pm 2$ °C bei einem Sollwert von 10 °C bei Umgebungstemperaturen von $\leq 25$ °C und $\leq 50$ % relativer Luftfeuchtigkeit
Gewicht	24 kg
Abmessungen (H x B x T)	$36 \times 42 \times 51$ cm
PC-Anschluss	Alle Funktionen und Parameter software-gesteuert über USB 1.1 Integrierter USB-Hub mit drei USB 1.1-Schnittstellen
E/A-Schnittstellen	4 digitale Eingänge, 4 programmierbare Relaisausgänge
Sicherheitsmerkmale	Leaksensor, automatische Platten-/Probenfläschchenerkennung, interne Überwachung aller mechanischen Operationen Überwachung der Funktion und Kühlleistung über 3 Temperatur- sensoren
Eingabe/Anzeige	LCD zur Anzeige von Systemparametern, Standby-Taste 3 LEDs (Power, Connected und Status) zur Statusangabe 4 Funktionstasten zum Betrieb während der Erstinstallation und zur Wartung

	WPS-3000TBRS
GLP	Unter Chromeleon: Automatische Gerätequalifikation (AutoQ <sup>™</sup> - Routinen) und Überwachung der System Wellness. Alle Systemparameter werden im Chromeleon Audit Trail protokolliert.
Medienberührte Teile	
Flusspfad Probe	Keramik, Titan, PEEK, MP35N
Flusspfad Waschflüssigkeit	Titan, PEEK, MP35N, PCTFE, Fused Silica, ETFE, ECTFE, PTFE, Glas
Leistungsaufnahme	100–120 V, 60 Hz; 200–240 V, 50 Hz; max. 320 VA
Emissionsschalldruckpegel	Typisch $< 65 \text{ dB}(A)$ in 1 m Abstand
Umgebungsbedingungen	Verwendungsbereich:InnenraumTemperaturbereich:10 °C bis 35 °CLuftfeuchtigkeit:80% relative Feuchte, nicht kondensierendÜberspannungskat.:IIVerschmutzungsgrad:2

# 9.5 WPS-3000TXRS

	WPS-3000TXRS
Probenkapazität	Abhängig von der Drehtellerkonfiguration:
Fläschchen	<ul> <li>216 x 0,3-mL-Fläschchen, 120 x 1,1-mL konische Fläschchen,</li> <li>216 x 1,2-mL-Fläschchen, 120 x 1,8- mL/2,0-mL-Fläschchen,</li> <li>66 x 4-mL-Fläschchen, 30 x 10- mL-Fläschchen bzw.</li> <li>120 x 0,5-mL oder 1,5-mL Eppendorf Fläschchen (offen, ohne Kappe)</li> <li>+ 15 x 10 mL Probenfläschchen</li> </ul>
Well Plates	3 x 96 bzw. 3 x 384 Mikrotiterplatten oder Deep-Well-Plates, 3 x 24 Deep-Well-Plates, 3 x 384 Low-Well-PCR-Plates + 15 x 10 mL Probenfläschchen
Injektionsmethode	Inline-Split-Loop
<b>Injektionsvolumen</b> Einstellbar (empfohlen) Optional	0,001 - 25 μL (0,2 – 25 μL) 0,1 – 100 μL
Genauigkeit Injektionsvolumen	Typisch ±0,5% bei 20 µL
Erforderliche Probenmenge (Minimum)	1 μL von 5 μL (250-μL-Fläschchen, konisch)
Präzision des Injektionsvolumens	${<}0,\!25$ % RSD bei 2 $\mu L$ (typisch ${<}0,\!15\%$ RSD), Koffein in Wasser
Linearität	Korr. koeff. >0,9999, RSD <0.5% bei 1-20 µL, Koffein in Wasser
Waschen der Nadel	Aktives Waschen der Probennadel von außen
Verschleppung	<0,004% für Koffein mit externem Waschen bei 20 MPa
Injektionszykluszeit	<20 s für 5 μL
Probenkühlung	4 bis 45 °C oder 22 °C unterhalb der Umgebungstemperatur
Genauigkeit Probentemperatur	$\pm 2$ °C bei einem Sollwert von 10 °C bei Umgebungstemperaturen von $\leq 25$ °C und $\leq 50$ % relativer Luftfeuchtigkeit
Gewicht	24 kg
Abmessungen (H x B x T)	$36 \times 42 \times 51$ cm
PC-Anschluss	Alle Funktionen und Parameter software-gesteuert über USB 1.1 Integrierter USB-Hub mit drei USB 1.1-Schnittstellen
E/A-Schnittstellen	4 digitale Eingänge, 4 programmierbare Relaisausgänge
Sicherheitsmerkmale	Leaksensor, automatische Platten-/Probenfläschchenerkennung, interne Überwachung aller mechanischen Operationen Überwachung der Funktion und Kühlleistung über 3 Temperatursensoren
Eingabe/Anzeige	LCD zur Anzeige von Systemparametern, Standby-Taste 3 LEDs (Power, Connected und Status) zur Statusangabe 4 Funktionstasten zum Betrieb während der Erstinstallation und zur Wartung

	WPS-3000TXRS
GLP	Unter Chromeleon: Automatische Gerätequalifikation (AutoQ <sup>™</sup> - Routinen) und Überwachung der System Wellness Alle Systemparameter werden im Chromeleon Audit Trail protokolliert.
Medienberührte Teile	
Flusspfad Probe	PEEK <sup>TM</sup> , Edelstahl, Vespel <sup>®</sup>
Flusspfad Waschflüssigkeit	PEEK, Edelstahl, PCTFE, Fused Silica, ETFE, ECTFE, PTFE, Glas
Leistungsaufnahme	100–120 V, 60 Hz; 200–240 V, 50 Hz; max. 320 VA
Emissionsschalldruckpegel	Typisch $< 65 \text{ dB}(A)$ in 1 m Abstand
Umgebungsbedingungen	Verwendungsbereich:InnenraumTemperaturbereich:10 °C bis 35 °CLuftfeuchtigkeit:80% relative Feuchte, nicht kondensierendÜberspannungskat.:IIVerschmutzungsgrad:2
# 10 Zubehör, Ersatzteile und Verbrauchsmaterialien

Zubehör, Ersatzteile und Verbrauchsmaterialien werden laufend dem neuesten technischen Standard angepasst. Eine Änderung der Bestellnummern ist deshalb nicht auszuschließen. Es ist jedoch sichergestellt, dass bei Bestellung der aufgeführten Bestellnummern stets voll kompatible Teile geliefert werden.

#### 10.1 Standardzubehör

Das folgende Standardzubehör ist im Lieferumfang enthalten (Änderungen vorbehalten). Einige der genannten Zubehörteile sind in den entsprechenden Ersatzteilkits enthalten. Informationen zum Inhalt der Kits finden Sie im Kapitel 10.3 ( $\rightarrow$  Seite 183).

Die Bestellnummer bezieht sich immer auf die jeweilige Verpackungseinheit. Wenn nicht anders angegeben, ist die Verpackungseinheit 1 Stück. Fragen hierzu beantwortet Ihnen gern die Thermo Fisher Scientific-Vertriebsgesellschaft für Dionex HPLC-Produkte.

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	1
Probenhalter (für 40 Probenfläschchen, A.D. < 12 mm) und Standard-Mikrotiterplatten (Höhe 12 - 24 mm)	6820.4070	3
Waschflasche WPS	Enthalten in 6820.4075	1
Sicherung, 2 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung, 4 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung 10 A, träge (6,3x 32 mm)	Enthalten in 6820.0026	1
Doppelmaulschlüssel (Größe 1/4 x 5/16)	6000.0051	1
Maulschlüssel (Größe 5,5 mm)	6146.1225	1
Maulschlüssel (Größe 3/16")	6146.1231	1
Sechskant-Stiftschlüssel (Größe 9/64")	6000.0053	1
Sechskant-Stiftschlüssel (Größe 2,5 mm)	6146.2625	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Probenfläschchen (1,8 mL), amber	Enthalten in 6000.0072	10
Bördelkappe mit Septum für 1,8-mL-Probenfläschchen	Enthalten in 6000.0071	10
Probenfläschchen (zylindrisch, 2 mL), amber für Schraubkappe	Enthalten in 6000.0060	10
Schraubkappe, weiß, für 2-mL-Probenfläschchen	Enthalten in 6000.0057	10

Bezeichnung	BestNr.	Menge im Zubehör
Silikondichtung, rot, für 2-mL-Probenfläschchen	Enthalten in 6000.0058	10
Probenfläschchen (zylindrisch, 10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Polypropylen-Probenfläschchen (250 µL)	Enthalten in 6820.0029	10
Deckel für 250-µL-Polypropylen-Probenfläschchen	Enthalten in 6820.0028	10
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000TBSL Analytical mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	1
Probenhalter (für 40 Probenfläschchen, A.D. < 12 mm) und Standard-Mikrotiterplatten (Höhe 12 - 24 mm)	6820.4070	3
Waschflasche WPS	Enthalten in 6820.4075	1
Sicherung, 2 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung, 4 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung 10 A, träge (6,3x 32 mm)	Enthalten in 6820.0026	1
Doppelmaulschlüssel (Größe 1/4 x 5/16)	6000.0051	1
Maulschlüssel (Größe 5,5 mm)	6146.1225	1
Maulschlüssel (Größe 3/16")	6146.1231	1
Sechskant-Stiftschlüssel (Größe 9/64")	6000.0053	1
Sechskant-Stiftschlüssel (Größe 2,5 mm)	6146.2625	1
Kapillare (PEEK, 0,18 mm x 1/16" ID x AD)	6827.5002	2
Fitting, fingertight, Rheflex (1/16", PEEK)	Enthalten in 6000.0012	4
Fittingschraube (1/16", 15 mm, PEEK)	6266.0024	3
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Probenfläschchen (1,8 mL), amber	Enthalten in 6000.0072	10
Bördelkappe mit Septum für 1,8-mL-Probenfläschchen	Enthalten in 6000.0071	10
Probenfläschchen (zylindrisch, 2 mL), amber für Schraubkappe	Enthalten in 6000.0060	10
Schraubkappe, weiß, für 2-mL-Probenfläschchen	Enthalten in 6000.0057	10
Silikondichtung, rot, für 2-mL-Probenfläschchen	Enthalten in 6000.0058	10

Bezeichnung	BestNr.	Menge im Zubehör
Probenfläschchen (zylindrisch, 10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Polypropylen-Probenfläschchen (250 µL)	Enthalten in 6820.0029	10
Deckel für 250-µL-Polypropylen-Probenfläschchen	Enthalten in 6820.0028	10
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000SL Semiprep, mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	1
Probenhalter (für 22 4-mL-Probenfläschchen)	6820.4084	3
Waschflasche WPS	Enthalten in 6820.4075	1
Sicherung, 2 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung, 4A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung 10A, träge (6,3x 32 mm)	Enthalten in 6820.0026	1
Doppelmaulschlüssel (Größe 1/4 x 5/16)	6000.0051	1
Maulschlüssel (Größe 5,5 mm)	6146.1225	1
Maulschlüssel (Größe 3/16")	6146.1231	1
Sechskant-Stiftschlüssel (Größe 9/64")	6000.0053	1
Sechskant-Stiftschlüssel (Größe 2,5 mm)	6146.2625	1
Probenfläschchen (4 mL, amber) für Schraubkappe	Enthalten in 6000.0074	10
Schraubkappe für 4-mL-Probenfläschchen	Enthalten in 6000.0073	10
Septum (Silikon/PTFE) für 4-mL-Probenfläschchen	Enthalten in 6000.0075	10
Probenfläschchen (zylindrisch, 10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000TBRS mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	1
Probenhalter (für 40 Probenfläschchen, A.D. < 12 mm) und Standard-Mikrotiterplatten (Höhe 12 - 24 mm)	6820.4070	3
Waschflasche WPS	Enthalten in 6820.4075	1
Sicherung, 2 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung, 4 A, träge (5 x 20 mm)	Enthalten in 6820.0026	2
Sicherung 10 A, träge (6,3x 32 mm)	Enthalten in 6820.0026	1
Doppelmaulschlüssel (Größe 1/4 x 5/16)	6000.0051	1
Maulschlüssel (Größe 5,5 mm)	6146.1225	1
Maulschlüssel (Größe 3/16")	6146.1231	1
Sechskant-Stiftschlüssel (Größe 9/64")	6000.0053	1
Sechskant-Stiftschlüssel (Größe 2,5 mm)	6146.2625	1
Kapillar-Kit (Viper) für UltiMate 3000 Bio-RS-Systeme (Single-Stack-Aufbau) mit 1 Kapillare (0,10 x 250 mm (ID x L), Viper, MP35N) z.B. für Verbindung von TCC-3000RS (Säulenauslass) zu DAD-3000RS, MWD-3000RS oder VWD-3400RS 1 Kapillare (0,10 x 350 mm (ID x L), Viper, MP35No) z.B. für Verbindung von WPS-3000TBRS zu TCC-3000RS (Säuleneinlass) 1 Kapillare (0,18 x 550 mm (ID x L), Viper, MP35No) z.B. für die Verbindung von RS-Pumpe zu WPS-3000TBRS	6841.2301	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Probenfläschchen (1,8 mL), amber	Enthalten in 6000.0072	10
Bördelkappe mit Septum für 1,8-mL-Probenfläschchen	Enthalten in 6000.0071	10
Probenfläschchen (zylindrisch, 2 mL), amber für Schraubkappe	Enthalten in 6000.0060	10
Schraubkappe, weiß, für 2-mL-Probenfläschchen	Enthalten in 6000.0057	10
Silikondichtung, rot, für 2-mL-Probenfläschchen	Enthalten in 6000.0058	10
Probenfläschchen (zylindrisch, 10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Polypropylen-Probenfläschchen (250 µL)	Enthalten in 6820.0029	10
Deckel für 250-µL-Polypropylen-Probenfläschchen	Enthalten in 6820.0028	10

Bezeichnung	BestNr.	Menge im Zubehör
Probenschleife, 100 µL (V=130 µL), MP35N, Viper	6841.2451	1
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 1 m)	6035.9035	1

## 10.2 Optionales Zubehör

Adapter	BestNr.
Adapter für 5-Positionen Halter zur Verwendung mit 2-mL-Probenfläschchen <i>Hinweis:</i> Bei Verwendung von 2-mL-Probenfläschchen im 5-Positionen Halter <i>muss</i> der Wert für <b>SampleHeightOffset_T</b> angepasst werden ( $\rightarrow$ Seite 92). Andernfalls wird die Probennadel beschädigt.	6820.4092
Adapter für Niederhalter Der Adapter wird für Deep-Well-Plates mit 24 Kavitäten benötigt. Bei Verwendung von 10-mL-Probenfläschchen ohne Kappe wird die Installation des Adapters ebenfalls empfohlen.	6820.2402
<i>Hinweis</i> : Wird der Adapter installiert, <i>muss</i> der automatische Tray-Test ausgeschaltet werden ( $\rightarrow$ Seite 94).	
Adapter für Low-Well-PCR-Plates (8 - 12 mm Plate-Höhe)	6820.4088

Injektionsvolumen-Kits		BestNr.
WPS-3000SL Analytical WPS-3000RS	250-μL-Kit, Viper, geeignet für Applikationen mit einem Druck <103 MPa (15000 psi) Das Kit enthält: Probenschleife (250 μL), SST, Viper Pufferschleife (>250 μL), SST, Viper Probennadel, SST, Viper-kompatibel Sicherungsschraube Probennadel Spritze (250 μL)	6822.2442
	<ul> <li>500-μL-Kit, Viper, geeignet für Applikationen mit einem Druck &lt;103 MPa (15000 psi)</li> <li>Das Kit enthält:</li> <li>1 Probenschleife (500 μL), SST, Viper</li> <li>1 Pufferschleife (&gt;250 μL), SST, Viper</li> <li>1 Probennadel, SST, Viper-kompatibel</li> <li>1 Sicherungsschraube Probennadel</li> <li>1 Spritze (500 μL)</li> </ul>	6822.2443
WPS-3000TBRS	<ul> <li>250-μL-Kit, Viper, geeignet für Applikationen mit einem Druck &lt;103 MPa (15000 psi)</li> <li>Das Kit enthält:</li> <li>1 Probenschleife (250 μL), MP35N, Viper</li> <li>1 Pufferschleife (&gt;250 μL), MP35N, Viper</li> <li>1 Probennadel, Keramik, Viper-kompatibel</li> <li>1 Sicherungsschraube Probennadel</li> <li>1 Spritze (250 μL)</li> <li>1 Nadelsitzkapillare (3,1 μL), MP35N, Viper</li> </ul>	6841.2442

Injektionsvolumen-Kits		BestNr.
WPS-3000TBRS (Fortsetzung)	<ul> <li>500-μL-Kit, Viper, geeignet für Applikationen mit einem Druck &lt;103 MPa (15000 psi)</li> <li>Das Kit enthält:</li> <li>1 Probenschleife (500 μL), MP35N, Viper</li> <li>1 Pufferschleife (&gt;250 μL), MP35N, Viper</li> <li>1 Probennadel, Keramik, Viper-kompatibel</li> <li>1 Sicherungsschraube Probennadel</li> <li>1 Spritze (500 μL)</li> <li>1 Nadelsitzkapillare (3,1 μL), MP35N, Viper</li> </ul>	6841.2443
WPS-3000SL Semiprep	<ul> <li>1000-μL -Kit für WPS-3000SL Semiprep</li> <li>Das Kit enthält:</li> <li>Probenschleife (1000 μL) und Spritze (1000 μL)</li> <li>Die Komponenten in diesem Kit sind für den WPS-3000SL Semiprep gedacht. Wenn Sie diese</li> <li>Komponenten in einem WPS-3000 Analytical</li> <li>verwenden möchten, ist ein Upgrade auf die</li> <li>semipräparative Autosamplerkonfiguration erforderlich.</li> <li>Wenden Sie sich in diesem Fall an den Thermo Fisher</li> <li>Scientific-Kundendienst für Dionex HPLC-Produkte.</li> </ul>	6822.2436

Probenhalter für	BestNr.
10 x 10-mL-Probenfläschchen und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4086
40 x 1,1-mL-Probenfläschchen ( <i>konisch</i> ) und/oder Standard-Mikrotiterplatten (Platten-Höhe 12 - 24 mm)	6820.4087
40 x 1,5-mL-Probenfläschchen ( <i>Eppendorf</i> ) Die Fläschchen müssen offen sein.	6820.4094
40 x 0,5-mL-Probenfläschchen ( <i>Eppendorf</i> ) Die Fläschchen müssen offen sein.	6820.4096
72 x 1,2-mL-Probenfläschchen ( <i>zylindrisch</i> ) und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4090
72 x 0,3-mL-Probenfläschchen ( <i>zylindrisch</i> ) und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4091
72 x 0,3 mL Mikrodialyse-Probenfläschchen	6820.4097
Deep-Well-Plate (Plate-Höhe: 34 - 46 mm)	6820.4079
Deep-Well-Plate (Plate-Höhe: 30 - 36 mm)	6820.4083
Deep-Well-Plate (Plate-Höhe: 20 - 32 mm)	6820.4089

Probenschleifen		BestNr.
WPS-3000SL Analytical, WPS-3000RS, WPS-3000TXRS	25 μL (V=40 μL), SST, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6820.2452
	100 μL (V=130 μL), SST, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6820.2451
	250 μL (V=344 μL), SST, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6820.2453
	500 μL (V=667 μL), SST, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6820.2454
WPS-3000TBRS	25 μL (V=40 μL), MP35N, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6841.2452
	100 μL (V=130 μL), MP35N, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6841.2451
	250 μL (V=344 μL), MP35N, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6841.2453
	500 $\mu$ L (V=667 $\mu$ L), MP35N, Viper Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6841.2454

Probenschleifen		BestNr.
WPS-3000SL Semiprep	1000 $\mu$ L (V=1618 $\mu$ L), SST Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. Diese Probenschleife ist für den WPS-3000SL Semiprep gedacht. Wenn Sie diese semipräparative Probenschleife in einem WPS-3000SL Analytical verwenden möchten, ist ein Upgrade auf die semipräparative Autosamplerkonfiguration erforderlich. Wenden Sie sich in diesem Fall an den Thermo Fisher Scientific-Kundendienst für Dionex HPLC-Produkte.	6820.2429

Spritzen	BestNr.
25 μL	6822.0001
250 μL	6822.0003
500 µL	6822.0004
1000 μL	6822.0005

Kapillar-Kits	BestNr.
Kit für UltiMate 3000 RSLC-Systeme (Single-Stack-Aufbau) mit LPG-3400RS oder DGP-3600RS Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,13 x 250 mm, 0,13 x 350 mm und 0,18 x 450 mm ID x L).	6040.2301
Kit für UltiMate 3000 RSLC-Systeme (Single-Stack-Aufbau) mit HPG-3200RS oder HPG-3400RS Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,13 x 250 mm, 0,13 x 350 mm und 0,18 x 550 mm ID x L).	6040.2308
Kit für UltiMate 3000 Bio-RS-Systeme (Single-Stack-Aufbau) mit UltiMate 3000 RS-Pumpe Das Kit enthält 3 Viper-Kapillaren, MP35N (je 1 x 0,10 x 250 mm, 0,10 x 350 mm und 0,18 x 550 mm (ID x L)).	6841.2301
Kit für UltiMate 3000 Standard-Systeme (Single-Stack-Aufbau) mit ISO-31000SD, LPG-3400SD oder DGP-3600SD Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,18 x 250 mm, 0,18 x 350 mm und 0,18 x 450 mm ID x L).	6040.2302
Kit für UltiMate 3000 Standard-Systeme (Single-Stack-Aufbau) mit HPG-3200SD oder HPG-3400SD Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,18 x 250 mm, 0,18 x 350 mm und 0,18 x 550 mm ID x L).	6040.2309
Kit für LPG-3400XRS, mit 2 Viper-Kapillaren, SST (je 1x 0,1 x 350 mm und 0,13 x 550 mm (ID x L)) 1 Viper-Kapillare, PEEK (0.065 x 250 mm (ID x L))	6043.2301

Sonstiges	BestNr.
Frontabdeckung, transparent Verschließt den Probenraum zum Schutz gegen Staub oder andere Partikel.	6820.1427
Signalkabel (6-pol. Mini-DIN) zum Anschluss externer Geräte an die Digital I/O-Buchse	6000.1004
Systemdrainage-Kit für UltiMate 3000 Systeme Das Kit enthält alle erforderlichen Komponenten für den Anschluss der System- drainage sowie eine detaillierte Installationsanleitung.	6040.0005
Upgrade-Kit WPS-3000SL Semiprep zur Umrüstung von WPS-3000SL Analytical auf WPS-3000SL Semiprep. Das Kit enthält alle für die Umrüstung erforderlichen Teile sowie eine 2500-µL-Proben- schleife. Das Upgrade wird vom Thermo Fisher Scientific-Kundendienst installiert.	6822.2450

#### **10.3 Ersatzteile und Verbrauchsmaterialien**

Die Bestellnummer bezieht sich immer auf die jeweilige Verpackungseinheit. Wenn nicht anders angegeben, ist die Verpackungseinheit 1 Stück.

Fragen hierzu beantwortet Ihnen gern die Thermo Fisher Scientific-Vertriebsgesellschaft für Dionex HPLC-Produkte.

Bezeichnung	BestNr.
Adapter für 5-Positionen Halter für die Verwendung von 2-mL-Probenfläschchen <i>Hinweis:</i> Bei Verwendung von 2-mL-Probenfläschchen im 5-Positionen Halter <i>muss</i> der Wert für <b>SampleHeightOffset_T</b> angepasst werden ( $\rightarrow$ Seite 92). Andernfalls wird die Probennadel beschädigt.	6820.4092
Adapter für Niederhalter Der Adapter wird für Deep-Well-Plates mit 24 Kavitäten benötigt. Bei Verwendung von 10-mL-Probenfläschchen ohne Kappe wird die Installation des Adapters ebenfalls empfohlen. <i>Hinweis</i> : Wird der Adapter installiert, <i>muss</i> der automatische Tray-Test ausgeschaltet werden ( $\rightarrow$ Seite 94).	6820.2402
Adapter für Low-Well-PCR-Plates (8 – 12 mm Plate-Höhe)	6820.4088
Diagnosetoolkit Das Kit wird zur Durchführung der Autosampler-Diagnosetests unter Chromeleon benötigt. Es enthält ein Verschlussfitting und eine Abfallkapillare.	6822.0030
Doppelmaulschlüssel (1/4 x 5/16)	6000.0051
Fitting (RH fingertight, 1/16", PEEK, einteilig), 10 Stück	6000.0012
Fitting- und Schneidringkit (RH) $\rightarrow$ Schneidring- und Fittingkit (RH)	
Fitting- und Schneidringkit (SR/FS-7) $\rightarrow$ Schneidring- und Fittingkit (SR/FS-7)	
Fittingschraube (PEEK, 1/16", 15 mm)	6266.0024
Injektionsvolumen-Kit (250 μL), Viper, für WPS-3000SL Analytical und WPS-3000RS geeignet für Applikationen mit einem Druck <103 MPa (15000 psi) Das Kit enthält: 1 Probenschleife (250 μL), SST, Viper 1 Pufferschleife (>250 μL), SST, Viper 1 Probennadel, SST, Viper-kompatibel, und 1 Sicherungsschraube Probennadel 1 Spritze (250 μL)	6822.2442
Injektionsvolumen-Kit (250 μL), Viper, für WPS-3000TBRS geeignet für Applikationen mit einem Druck <103 MPa (15000 psi) Das Kit enthält: 1 Probenschleife (250 μL), MP35N, Viper 1 Pufferschleife (>250 μL), MP35N, Viper 1 Probennadel, Keramik, Viper-kompatibel und 1 Sicherungsschraube Probennadel 1 Spritze (250 μL) 1 Nadelsitzkapillare (3,1 μL), MP35N, Viper	6841.2442

Bezeichnung	BestNr.
Injektionsvolumen-Kit (500 μL), Viper, für WPS-3000SL Analytical und WPS-3000RS geeignet für Applikationen mit einem Druck <103 MPa (15000 psi) Das Kit enthält: 1 Probenschleife (500 μL), SST, Viper 1 Pufferschleife (>250 μL), SST, Viper 1 Probennadel, SST, Viper-kompatibel, und 1 Sicherungsschraube Probennadel 1 Spritze (500 μL)	6822.2443
Injektionsvolumen-Kit (500 μL), Viper, für WPS-3000TBRS geeignet für Applikationen mit einem Druck <103 MPa (15000 psi) Das Kit enthält: 1 Probenschleife (500 μL), MP35N, Viper 1 Pufferschleife (>250 μL), MP35N, Viper 1 Probennadel, Keramik, Viper-kompatibel und 1 Sicherungsschraube Probennadel 1 Spritze (500 μL) 1 Nadelsitzkapillare (3,1 μL), MP35N, Viper	6841.2443
Injektionsvolumen-Kit (1000 μL), WPS-3000SL Semiprep Das Kit enthält: Probenschleife (1000 μL) und Spritze (1000 μL) Die Komponenten in diesem Kit sind für den WPS-3000SL Semiprep gedacht. Wenn Sie diese Komponenten in einem WPS-3000SL Analytical verwenden möchten, ist ein Upgrade auf die semipräparative Autosamplerkonfiguration erforderlich. Wenden Sie sich in diesem Fall an den Thermo Fisher Scientific- Kundendienst.	6822.2436
Injektor-Waste-Kapillare, Viper, für WPS-3000SL, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS	6820.2469
Kapillare, SST, Viper (0,18 x 450 mm ID x L) Verbindung Pumpe - WPS-3000SL Analytical	6040.2365
Kapillare, SST (0,25 x 1,58 mm ID x AD) Verbindung WPS-3000SL Analytical – TCC-3000	6820.2418
Kapillare, PEEK (90 μm x 250 mm ID x AD) z.B. für die Verbindung TCC-3000 (Säulenausgang) – Detektor in UltiMate 3000 Bio-RS-Systemen	6041.9025
Kapillare, PEEK (0,18 mm x 1/16" ID x AD), WPS-3000TBSL Analytical	6827.5002
Kapillare, MP35N, Viper (0,10 x 250 mm ID x L) z.B. für die Verbindung vom TCC-3000RS (Säulenauslass) zum DAD-3000RS, MWD-3000RS oder VWD-3400RS	6042.2330
Kapillare, MP35N, Viper (0,10 x 350 mm ID x L) z.B. für die Verbindung vom WPS-3000TBRS zum TCC-3000RS (Säuleneinlass)	6042.2340
Kapillare MP35N, Viper (0,18 x 550 mm (ID x L) z.B. für die Verbindung von RS-Pumpe zum WPS-3000TBRS	6042.2355
Kapillar-Kit für UltiMate 3000 RS-System mit LPG-3400RS oder DGP-3600RS Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,13 x 250 mm, 0,13 x 350 mm und 0,18 x 450 mm ID x L).	6040.2301

Bezeichnung	BestNr.
Kapillar-Kit für UltiMate 3000 RS-System mit HPG-3200RS oder HPG-3400RS Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,13 x 250 mm, 0,13 x 350 mm und 0,18 x 550 mm ID x L).	6040.2308
Kapillar-Kit für UltiMate 3000 Bio-RS-System mit UltiMate 3000 RS-Pumpe Das Kit enthält 3 Viper-Kapillaren, MP35N (je 1 x 0,10 x 250 mm, 0,10 x 350 mm und 0,18 x 550 mm (ID x L)).	6841.2301
Kapillar-Kit für UltiMate 3000 SD-System mit ISO-31000SD, LPG-3400SD oder DGP-3600SD Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,18 x 250 mm, 0,18 x 350 mm und 0,18 x 450 mm ID x L).	6040.2302
Kapillar-Kit für UltiMate 3000 SD-System mit HPG-3200SD oder HPG-3400SD Das Kit enthält 3 Viper-Kapillaren (SST; je 1x 0,18 x 250 mm, 0,18 x 350 mm und 0,18 x 550 mm ID x L).	6040.2309
Kapillarkit für LPG-3400XRS Das Kit enthält 2 Viper-Kapillaren, SST (je 1x 0,1 x 350 mm und 0,13 x 550 mm (ID x L)) 1 Viper-Kapillare, PEEK (0.065 x 250 mm (ID x L))	6043.2301
Maulschlüssel (Größe 5,5 mm)	6146.1225
Maulschlüssel (Größe 3/16")	6146.1231
Menüstift	6300.0100
Mikrotiterplatte (Standard-Mikrotiterplatte mit 96 Kavitäten)	6820.4100
Mikrotiterplatte (Deep-Well-Plate mit 96 Kavitäten)	6820.4101
Mikrotiterplatte (Standard-Mikrotiterplatte mit 384 Kavitäten)	6820.4110
Mikrotiterplatte (Deep-Well-Plate mit 384 Kavitäten)	Auf Anfrage
Nadelsitz, 2 Stück, für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS WPS-3000TBRS	6820.0047A 6841.0047
Nadelsitz-Servicekit mit Nadelsitz und Halteschraube für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS WPS-300TBRS WPS-3000SL Semiprep	6820.0038A 6841.0038 6820.0043
Nadelsitzkapillare, 3,1 μL, SST, Viper, für WPS-3000SL Analytical und WPS-3000RS <i>Hinweis</i> : Die Nadelsitzkapillare ist für Applikationen mit einem Druck <125 MPa	6820.2464
(18130 psi) geeignet.	
Nadelsitzkapillare, 3,1 µL, PEEK, Viper, für WPS-3000TBSL Analytical	6827.2408
Nadelsitzkapillare, 3,1 μL, MP35N, Viper, für WPS-3000TBRS <i>Hinweis</i> : Die Nadelsitzkapillare ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6841.2464
Nadelsitzkapillare, 1 µL, MP35N, Viper, für WPS-3000TBRS <i>Hinweis</i> : Die Nadelsitzkapillare ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6841.2472

Bezeichnung	BestNr.
Nadelsitzkapillare, 1 µL, SST, Viper, fürWPS-3000TXRS <i>Hinweis</i> : Die Nadelsitzkapillare ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6820.2472
Nadelsitzkapillare, 24 µL, SST, für WPS-3000SL Semiprep	6820.2409
Netzkabel, EU	6000.1000
Netzkabel, US	6000.1001
Netzkabel, UK	6000.1020
Netzkabel, Schweiz	6000.1030
Netzkabel, Italien	6000.1040
Netzkabel, Japan	6000.1050
Netzkabel, Australien, China	6000.1060
Netzkabel, Dänemark	6000.1070
Netzkabel, Indien/SA	6000.1090
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450
Pod (einschl. Rotor Seal für pH-Werte 0-14) für Injektionsventil imWPS-3000SL Analytical/Semiprep (Druck < 62 MPa (9000 psi))	6822.0051 6723.9013 6840.0013 6841.0013
Probenfläschchen (zvlindrisch 250 µL: Polypropylen: 1000 Stück)	6820.0029
Probenfläschchen (konisch. 1.1 mL. für Bördelkappe, 500 Stück)	6000.0077
Probenfläschchen (zvlindrisch, 1.2 mL, für Bördelkappe, 500 Stück)	6000.0062
Probenfläschchen (1.8 mL. braun, 100 Stück)	6000.0072
Probenfläschchen (2 mL; braun, für Schaubkappe; 500 Stück)	6000.0060
Probenfläschchen (4 mL; braun, für Schraubkappe; 100 Stück)	6000.0074
Probenfläschchen (zylindrisch, 10 mL), mit Bördelkappe und Septum (5 Stück)	6820.0023
Probenfläschchendeckel für 250-µL-Polypropylen-Fläschchen (1000 Stück)	6820.0028
Probenfläschchendeckel für 250-µL-Polypropylen-Fläschchen (100 Stück)	6820.0046
Probenfläschchendeckel für 1.1-mL-Fläschchen (Crimpkappe mit geschlitztem Silikon/PTFE-Septum; 100 Stück)	6000.0076
Probenfläschchendeckel für 1,2-mL-Fläschchen (Crimpkappe; 1000 Stück)	6000.0064
Probenfläschchendeckel mit Septum für 1,8-mL-Fläschchen (Crimpkappe; 100 Stück)	6000.0071

Bezeichnung	BestNr.
Probenfläschchendeckel (Schraubkappe, weiß) für 2-mL-Probenfläschchen (500 Stück)	6000.0057
Probenfläschchendeckel (Schraubkappe) für 4-mL-Probenfläschchen (100 Stück)	6000.0073
Probenfläschchenseptum (Silikon, rot) für 2-mL-Fläschchen (500 Stück)	6000.0058
Probenfläschchenseptum (Silikon/PTFE) für 4-mL-Probenfläschchen (500 Stück)	6000.0075
Probenfläschchenseptum (Silikon/PTFE, 8 mm, geschlitzt) für 1,2-mL-Probenfläschchen (500 Stück)	6000.0061
Probenhalter für 40 Probenfläschchen, A.D. < 12 mm und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4070
Probenhalter für 40 x 1,5-mL-Probenfläschchen ( <i>Eppendorf</i> ) Die Fläschchen müssen offen sein.	6820.4094
Probenhalter für 40 x 0,5-mL-Probenfläschchen ( <i>Eppendorf</i> ) Die Fläschchen müssen offen sein.	6820.4096
Probenhalter für 40 <i>konische</i> 1,1-mL-Probenfläschchen und/oder Standard- Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4087
Probenhalter für 72 zylindrische 1,2-mL-Probenfläschchen und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4090
Probenhalter für 72 x 0,3-mL-Probenfläschchen (zylindrisch) und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4091
Probenhalter für 72 x 0,3-mL-Mikrodialyse-Probenfläschen	6820.4097
Probenhalter für 10 x 10-mL-Probenfläschchen und/oder Standard- Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4086
Probenhalter für 22 x 4-mL-Probenfläschchen und/oder Standard-Mikrotiterplatten (Platten-Höhe: 12 - 24 mm)	6820.4084
Probenhalter für Deep-Well-Plate (Plate-Höhe: 30 - 36 mm)	6820.4083
Probenhalter für Deep-Well-Plate (Plate-Höhe: 34 - 46 mm)	6820.4079
Probenhalter für Deep-Well-Plate (Plate-Höhe: 20 - 32 mm)	6820.4089
Halter (5-Positionen) für 10-mL-Probenfläschchen Adapter für 5-Positionen-Halter für die Verwendung von 2-mL-Probenfläschchen	6820.4073 6820.4092
Probennadel, SST, Viper-kompatibel, für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS	6820.2432
Probennadel, Keramik, für WPS-3000TBRS	6841.2420
Probennadel, SST, für WPS-3000TBSL Analytical	6820.2403
Probennadel, SST, für WPS-3000SL Semiprep	6820.2419
Probennadel-Sicherungsschraube (Viper und Viper-kompatibel) für WPS-3000SL Analytical, WPS-3000SL Semiprep, WPS-3000RS, WPS-3000TBRS und WPS-3000TXRS	6820.2404

Bezeichnung	BestNr.
Probenschleife, Mikro, 25 μL (V=40 μL), SST, Viper für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6820.2452
Probenschleife, Mikro, 25 μL (V=40 μL), MP35N, Viper, für WPS-3000TBRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6841.2452
Probenschleife, analytisch, 100 μL (V=130 μL), SST, Viper für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6820.2451
Probenschleife, 100 µL (V=130 µL), MP35N, Viper, für WPS-3000TBRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <125 MPa (18130 psi) geeignet.	6841.2451
Probenschleife, analytisch, 100 µL (V=150 µL), PEEK, für WPS-3000TBSL Analytical Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen.	6820.2431
Probenschleife, 250 μL (V=344 μL), SST, Viper, für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6820.2453
Probenschleife, 250 µL (V=344 µL), MP35N, Viper, für WPS-3000TBRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6841.2453
Probenschleife, analytisch, 500 µL (V=667 µL), SST, Viper, für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6820.2454
Probenschleife, 500 μL (V=667 μL), MP35N, Viper für WPS-3000TBRS Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. <i>Hinweis</i> : Die Probenschleife ist für Applikationen mit einem Druck <103 MPa (15000 psi) geeignet.	6841.2454

Bezeichnung	BestNr.
Probenschleife, 1000 µL (V=1618 µL), SST, für WPS-3000SL Semiprep, Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. Diese Probenschleife ist für den WPS-3000SL Semiprep gedacht. Wenn Sie diese semipräparative Probenschleife in einem WPS-3000SL Analytical verwenden möchten, ist ein Upgrade auf die semipräparative Autosamplerkonfiguration erforderlich. Wenden Sie sich in diesem Fall an den Thermo Fisher Scientific- Kundendienst.	6820.2429
Probenschleife, 2500 µL (V=3300 µL), SST, für WPS-3000SL Semiprep Der Wert vor der Klammer gibt das zulässige Injektionsvolumen an. Der Wert in der Klammer ist das Gradientenverzögerungsvolumen. Diese Probenschleife ist für den WPS-3000SL Semiprep gedacht. Wenn Sie diese semipräparative Probenschleife in einem WPS-3000SL Analytical verwenden möchten, ist ein Upgrade auf die semipräparative Autosamplerkonfiguration erforderlich. Wenden Sie sich in diesem Fall an den Thermo Fisher Scientific- Kundendienst.	6820.2416
Pufferschleife, 100 µL, SST, Viper, für WPS-3000SL Analytical, WPS-3000RS und WPS-3000TXRS	6820.2466
Pufferschleife, 100 µL, MP35N, Viper, für WPS-3000TBRS	6841.2466
Pufferschleife, >250 µL, MP35N, Viper, für WPS-3000TBRS	6841.2468
Pufferschleife, >250 µL, SST, Viper, für WPS-3000SL Semiprep	6820.2468
Reinigungsstäbchen (10 Stück)	6040.0006
Rotor Seal (geeignet für pH-Werte 0 - 14) für Injektionsventil im WPS-3000SL Analytical/Semiprep ,WPS-3000RS und WPS-3000TBRS	6840.0012
Rotor Seal (geeignet für pH-Werte 0-10) für Injektionsventil im WPS-3000TXRS <i>Empfehlung</i> : Nach einem Tausch des Rotors Seals wird unter Umständen nicht immer die volle Lebensdauer des Ventils erreicht. Thermo Fisher Scientific empfiehlt daher, den Ventil-Pod zu tauschen ( $\rightarrow$ Seite 154).	6843.0012
Rotor Seal und Stator Face Seal kit (geeignet für pH-Werte 0-14) für Injektionsventil im WPS-3000TBSL Analytical	6722.9014
Schlauch zwischen Waschflasche und Spritzenventil	6820.0011
Sechskant-Stiftschlüssel (9/64")	6000.0053
Sechskant-Stiftschlüssel (2,5 mm)	6146.2625
Signalkabel (6-pol. Mini-DIN) zum Anschluss externer Geräte an die Digital I/O-Buchse	6000.1004
Sicherungen, Kit mit: Sicherung, 2 A, träge, 5x20 mm (10 St.) Sicherung, 4 A, träge, 5x20 mm (10 St.) Sicherung, 10 A, träge, 250V, 6,3 x 32 mm, 250 V (5 St.)	6820.0026
Sicherungsschraube Probennadel $\rightarrow$ Probennadel-Sicherungsschraube	
Spritze (25 µL)	6822.0001
Spritze (100 µL)	6822.0002
Spritze (250 µL)	6822.0003

Bezeichnung	BestNr.
Spritze (500 µL)	6822.0004
Spritze (1000 µL)	6822.0005
Spritze (2500 µL)	6822.0006
Spritzendichtung (zum Einbau in das Spritzenventil; 5 Dichtungen)	6822.0009
Stator für Injektionsventil im WPS-3000SL Analytical/Semiprep und WPS-3000RS WPS-3000TXRS <sup>1</sup> WPS-3000TBRS WPS-3000TBSL Analytical <sup>1</sup> <i>Empfehlung</i> : Nach einem Tausch des Stators wird unter Umständen nicht immer die volle Lebensdauer des Ventils erreicht. Thermo Fisher Scientific empfiehlt daher, den Ventil-Pod zu tauschen ( $\rightarrow$ Seite 154).	6840.0011 6840.0011 6841.0011 6723.9011
Systemdrainage-Kit für UltiMate 3000 Systeme Das Kit enthält alle erforderlichen Komponenten für den Anschluss der Systemdrainage sowie eine detaillierte Installationsanleitung.	6040.0005
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 5 m)	6911.0002
USB-Kabel, Typ A auf Typ B, High-speed USB 2.0 (Kabellänge: 1 m)	6035.9035
Waschflasche (2 Stück), einschließlich Halteklammer und Schlauch	6820.4075

# 11 Anhang - Anschlussbelegung Digital I/O

Über die Digital I/O-Buchsen stehen vier digitale Eingänge und vier Relaisausgänge zur Verfügung, die zur Übertragung von digitalen Signalen mit externen Geräten verwendet werden können. Die Nummerierung der Buchsen auf der Geräterückseite entspricht der Nummerierung der entsprechenden Relais und digitalen Eingänge in Chromeleon.



COM ist der gemeinsame Kontakt für NO und NC. Ist das Relais nicht aktiviert oder ist der Autosampler ausgeschaltet, besteht eine Verbindung zwischen COM und NC. Ist das Relais aktiviert, besteht eine Verbindung zwischen COM und NO.

\* NO = Normally Open (Arbeitskontakt) NC = Normally Closed (Ruhekontakt)

Abb. 89: Digital I/O (Ansicht: Geräterückseite)

	Beschreibung
Digital Input $\leftrightarrow$ Digital GND	0 bis +5V
$Relay\_Com \leftrightarrow Relay\_NO (bzw. Relay\_NC)$	Schaltspannung: 100V DC; Schaltstrom: 0,25 A Dauergrenzstrom: 0,5 A; Schaltleistung 3W Durchgangswiderstand: max. 200mOhm

Abb. 90: Digital I/O

Verwenden Sie zum Anschluss von externen Geräten das 6-adrige Mini-DIN-Signalkabel (Best.-Nr. 6000.1004). Informationen zur Funktion der Pins sowie zur Farbe und zum Label der Kabeladern, die mit dem jeweiligen Pin verbunden sind, finden Sie in der unten stehenden Tabelle.

**i** Hinweis: Wenn Sie eine Pumpe des Typs LPG-3400XRS anschließen möchten, finden Sie alle erforderlichen Information in der *Bedienungsanleitung* zur Pumpe.

Pin	Signalname	Signalpegel	Farbe Ader	Label Ader
1	nicht belegt		pink	Analog High
2	Relay_NC	max. 100V <sub>DC</sub> /0,25A/3W	grau	Analog Low
3	Digital_GND	0V	grün	Digital GND
5	Digital Input	0 bis +5V	gelb	Digital Input
6	Relay_NO	max. 100V <sub>DC</sub> /0,25A/3W	braun	Digital Output
8	Relay_COM	max. 100V <sub>DC</sub> /0,25A/3W	weiß	Dig./Out (GND)
			schwarz	Shield

Abb. 91: Anschlussbelegung (6-adrige Mini-DIN Buchse und Kabel)

# 12 Index

#### A

Anschluss	
Chromeleon-Rechner	35
Digital I/O	36, 191
Fluidik	26
Netzkabel	
Signalkabel	
Spritzenventil	53
USB	35
Waschfllüssigkeit	50
Anzeige	20
Äquilibrieren	60
Äquilibrieren (SmartStartup Wizard)	60
Äquilibrierprogramm	60
Auspacken	32
Außerbetriebnahme	98
Ausstattungsvarianten	14
Autosampler Drainage	54

#### B

Batch72Bedienelemente20Bedienungsanleitung1Beleuchtung25Betrieb63Anzeige20Bedienelemente20Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sich erk siteme fürstlande74	Back	76
Bedienelemente20Bedienungsanleitung1Beleuchtung25Betrieb63Anzeige20Bedienelemente20Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sich erk siteme fürstlande74	Batch	72
Bedienungsanleitung1Beleuchtung25Betrieb63Anzeige20Bedienelemente20Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Siehertenisteren63	Bedienelemente	
Beleuchtung25Betrieb63Anzeige20Bedienelemente20Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sich erk siteme fürstlande	Bedienungsanleitung	1
Betrieb    63      Anzeige    20      Bedienelemente    20      Chromeleon    67      Einschalten    63      Einstellungen    81      Empfehlungen    87      Funktionstasten    73      Gerätedisplay    73      Menü    75	Beleuchtung	25
Anzeige20Bedienelemente20Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sich erheitere fürstlande74	Betrieb	63
Bedienelemente20Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sich erheitere fürstlande74	Anzeige	20
Chromeleon67Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sichertheiterre74	Bedienelemente	
Einschalten63Einstellungen81Empfehlungen87Funktionstasten73Gerätedisplay73Menü75Sichertheiterre74	Chromeleon	67
Einstellungen	Einschalten	63
Empfehlungen	Einstellungen	
Funktionstasten   73     Gerätedisplay   73     Menü   75	Empfehlungen	
Gerätedisplay	Funktionstasten	73
Menü	Gerätedisplay	73
Ciahanhaitan a Quahanan (	Menü	75
Sicherneitsmabhanmen4	Sicherheitsmaßnahmen	4
Bypass-Modus13	Bypass-Modus	13

#### С

Cancel	77
Chromeleon	
Autosampler installieren	
Autosampler konfigurieren	40
Installationsprogramm	
Lizenz	
Predictive Performance	
Program Wizard	71
Programm erstellen	71
Server Configuration Program	
Server Monitor	

Spezielle Funktionen	. 88
Steuerung	. 67
Steuerung automatisch	. 71
Steuerung direkt	. 68
USB-Treiber	. 37
Verbinden	. 67
Chromeleon-Dialoge	
Error Levels	. 45
General	. 40
Inputs	. 44
Options	. 43
Relays	. 44
Segments / Pump Link	. 41
Sharing	. 41
Commands (Dialogfenster)	. 68
Configuration (Menü)	. 80
Control (Menü)	. 78
Control Panel	. 69
Coolant Temperature	. 79
Cooling Power	. 79
Cooling Pump Power	. 79

#### D

DCMSLink	
Deep-Well-Plates	
Device View	
Diagnose	
Diagnose-Meldungen	
Diagnostics (Menü)	79
Dialogfenster Commands	
Digital I/O	
Dispense Destination	
Display & Soft Keys	
Drainage	
Anschluss	54
Autosampler	54
Drehteller	
Konfiguration ändern	
Tray Shake	
Tray-Test	
2	

#### Е

Einschalten	63
Elektrischer Anschluss	
Entgasen	
Entlüften	
Error Levels (Dialogfenster)	
Ersatzteile	
Execute	77

#### F

Fan Power	79
Fehlermeldungen	
Fehlersuche	
Diagnose-Meldungen	
Fehlermeldungen	
Störungen	
Firmware aktualisieren	
Firmware-Download	
Fluidische Anschlüsse	
Funktionsprinzip	
Funktionstaste	73
Back	
Cancel	77
Clear	
Execute	77
Menu	74
Next	77, 104
Ok	77
Pfeil nach oben	
Pfeil nach rechts	
Pfeil nach unten	
Prev	77, 104
Prime	74
Rotate	74
Select	76
Wash	74

#### G

#### H

Heat Sink Temperature79
-------------------------

I

Inbetriebnahme	47
5-Positionen Halter	59
Allgemein	47

Anschluss Waschflüssigkeit	50
Äquilibrieren	60
Entlüften	55
Kapillaren anschließen	
Probenhalter	57
Spritze spülen	55
Spritzenventil	53
Inject Valve	
Inject Wash	80
Injektion (Synchronisation)	
Injektionsbefehl	
Injektionsparameter	84
Injektionsventil	
Injizieren	
Inputs (Dialogfenster)	44
Installation	
in DCMSLink	
Probenhalter	
Probennadel	
Probenschleife	
Pufferschleife	
Spritzenwechsel	
Standort	
unter Chromeleon	
Verbinden	

#### K

Kapillaren (Viper)	48
Kapillaren anschließen	48
Kommandos (Display)	73

#### L

Leak Sensor Mode	80, 86
Leak Test	
Leakerkennung	86
Leaksensor	26, 86, 120
Lösungsmittel	

#### Μ

Main (Menü)	
Maintenance (Menü)	
Meldungen	
Menü	
Aufbau	
Configuration	
Control	
Diagnostics	
Main	
Maintenance	
Übersicht	
Mikrotiterplatte	
Einsetzen	
Empfehlung	
Erkennung	
Module Address	

#### Ν

91
139
144
79
22
78

#### 0

Ok	77
Operational Qualification	94
Optimieren	161
Options (Dialogfenster)	43

#### Р

9
4
0
9
5
9
4
4
5
4
3
7
7
5
5
0
5
7
5
1
3
7
6
1
0
1
6
6
9
6
9
9
1
2
4
· •
1

#### R

Real Temperature	
Relative Probenposition	
Relays (Dialogfenster)	
Reset to Factory Defaults	80
RMA-Nummer	
Rotor Seal	

### S

Sample Cooling	
SampleHeight	
SampleHeightOffset	
Segments / Pump Link (Dialogfenster)	
Select	
Self Test	
Server Configuration (Programm)	
Server Configuration Program	39
Server Monitor	39
Service	
Allgemein	119
Firmware aktualisieren	159
Iniektionsventil	148
Leaksensor	120
Nadelsitz wechseln	139
Nadelsitzkanillare wechseln	144
Probennadel waschen	130
Probennadel wechseln	130
Probenschleife	125
Pufferschleife	124
Sicherheitsmaßnahmen	110
Sicherungswechsel	119
Spritzenwechsel	138
Spinzenweenser	121
Shutdown Programm	00
Sindaowii-1 logramm	
Sicherheitemeßnehmen	
Sichemin aaachlitten	4, 119
Sicherungsschnlitten	
Sicherungswechser	138
Signaikadei	
SmartShuldown	87, 99
	00, 87
Solikeys	
	103
Spritze	17, 121
Spritze spulen	
Spritzenventil	
Spulen (Spritze)	
Standby-Programm	
Stator	154
Stator Face Seal	155
Statusanzeige	
Display	
LEDs	
Status-LEDs	103

#### Steuerung

Automatisch	71
Chromeleon	67
Direkt	
Störungen	
Symbole	3
Synchronisation (Injektion)	
Syringe	
Systemdrainage	54

#### Т

Technische Daten	
Tray	23, 57, 80
Tray Shake	
Tray-Test	
5	

#### U

Udp-Kommando	96
USB	
USB-Anschluss	
USB-Konfigurationsdatei	

#### V

Verbindung (Autosampler-Pumpe)	
Verbrauchsmaterialien	
Version Info	79

# Verwendungszweck.8Viper-Kapillaren.48Vorbereitung.475-Positionen Halter59Allgemein47Anschluss Waschflüssigkeit50Äquilibrieren.60Entlüften55Kapillaren anschließen48Spritze spülen55Spritzenventil53

#### W

Wartung	
Wartungsintervalle	
Waschen (Probennadel)	
Waschflasche	
Waschflüssigkeit	
Wash	
Wellness	
Well-Plate	

#### Z

Zubehör	
Optional	178
Standardzubehör	173