



# Thermo Scientific Dionex UltiMate 3000-Serie

## Autosampler WPS-3000PL (RS) und WPS-3000FC

## Bedienungsanleitung (Originalbedienungsanleitung)



Version: 1.6 Datum: September 2013

© 2013 Thermo Fisher Scientific Inc. Dok-Nr. 4828.2001

# CE

# EG KONFORMITÄTSERKLÄRUNG

(Original-Konformitätserklärung)

Geräteart: Thermo Scientific Dionex UltiMate 3000 - Autosampler

Typenbezeichnung: WPS-3000TBPL WPS-3000PL RS und WPS-3000TPL RS WPS-3000TFC und WPS-3000TBFC

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 f
ür elektrische Mess-, Steuer-, Regel- und Laborger
äte 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 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	Sicherheit	3
	1.2	2.1 Symbole am Gerät und in der Bedienungsanleitung	3
	1.2	2.2 Sicherheitsmaßnahmen	4
	1.3	Verwendungszweck des Gerätes	8
า	ŕn	h out light	11
2	U	DerDIICK	•••• 11
	2.1	Kurzbeschreibung	11
	2.2	Funktionsprinzip	12
	2.3	Autosamplerkonfigurationen	13
	2.3	3.1 Übersicht	13
	2.3	3.2 Biokompatible Autosampler	14
	2.3	3.3 Autosampler mit Probenthermostatisierung	15
	2.3	3.4 Autosampler zur Fraktionssammlung	15
	2.3	3.5 Kombinationsmöglichkeiten von Spritzen und Schleifen	16
	2.4	Innenansicht	17
	2.4	4.1 WPS-3000PL RS und WPS-3000TBPL	17
	2.4	4.2 WPS-3000TFC und WPS-3000TBFC	19
	2.5	Gerätevorderseite	21
	2.6	Geräterückseite	22
	2.6	6.1 Netzschalter	23
	2.6	6.2 Sicherungsschlitten	23
	2.6	6.3 USB-Anschluss	23
	2.6	6.4 Digital I/O	23
	2.7	Drehteller, Probenhalter, Probengefäße und Mikrotiterplatten	24
	2.8	Probenraumbeleuchtung	27
	2.9	Fluidische Anschlüsse	27
	2.10	Leaksensor	27
	2.11	Injektionsventil	28
	2.12	Steuerung über Chromeleon	29
	2.13	Wellness und Predictive Performance	30
3	In	ostallation	31
5	2 1		
	3.1	Anforderungen an den Standort	31
	3.2	Auspacken	31
	3.3	Position des Autosamplers im UltiMate 3000-System	
	3.4	Verbinden des Autosamplers	34
	3.4	4.1 Allgemeine Informationen	34
	3.4	4.2 Anschluss des USB-Kabels	34
	3.4	4.3 Anschluss des Netzkabels	35
	3.4	4.4 Anschluss des Signalkabels (Digital I/O)	35

	3.5 Ein	richten des Autosamplers in Chromeleon	
	3.5.1	Laden des USB-Treibers für den Autosampler	
	3.5.2	Installieren des Autosamplers	40
	3.5.3	Konfigurieren des Autosamplers	41
	3.6 Ein	richten des Autosamplers in DCMSLink	48
4	Vorbe	reitung für den Betrieb (Inbetriebnahme)	49
	4.1 Übe	ersicht	49
	4.2 Hin	weise zum Anschluss von Kapillaren	51
	4.3 Ans	chließen der Waschflüssigkeit	52
	4.3.1	Anschluss der Waschflüssigkeit bei Online-Entgasung	52
	4.3.2	Anschluss der Waschflasche	54
	4.4 Ans	chlüsse am Spritzenventil	55
	4.5 Ans	chließen der Systemdrainage	56
	4.6 Ent	fernen von Luftblasen aus der Spritze (Prime)	56
	4.7 Ein	setzen und Bestücken eines Probenhalters	
	4.8 Bes	tücken der 5-Positionen-Halter	61
	4.9 Aqı	illibrieren des Systems	62
5	Betrie	b und Wartung	63
	5.1 Ein	schalten des Autosamplers	63
	5.2 Stat	usanzeige	64
	5.3 Pro	benposition	65
	5.4 Stee	aerung über Chromeleon	67
	5.4.1	Verbinden mit Chromeleon	67
	5.4.2	Direkte Steuerung	68
	5.4.3	Automatische Steuerung	71
	5.5 Fun	ktionstasten und Menüs am Gerätedisplay	74
	5.5.1	Einblenden der Funktionstasten	74
	5.5.2	Autosampler-Menüs	7/6
	5.6  Eins	stellungen für den Betrieb	83
	5.0.1	Festiegen der injektionsmetnode	83
	5.6.2	Einschalten der Drobenthermostetisierung	00
	5.0.5	Ändern der Drehtellerkonfiguration	00 97
	5.6.5	Fastlagen der Injektionsparameter und Starten der Injektion	07
	5.6.5	Finschalten der Leakerkennung	00
	5.6.7	Annassen von Helligkeit und Kontrast der Displayanzeige	9 89
	5.6.8	Empfehlungen für den Betrieb	90
	5.7 Sne	zielle Funktionen in Chromeleon	90
	571	Aktive Überwachung von Verschleißteilen (Predictive Performance)	91
	5.7.2	Definieren der Nadelhöhe und Vorstecher-Einstechtiefe	
	5.7.3	Bewegen des Drehtellers vor dem Aufziehen der Probe (Trav Shake)	94
	5.7.4	Operational Qualification und Performance Qualification	95
	5.7.5	Automatischer Tray-Test	95
	5.7.6	Automatische Probengefäßerkennung	95

5	5.7.7 Benutzerdefinierte Programme	
5.8	Außerbetriebnahme des Autosamplers	
5.9	Wartung und Wartungsintervalle	100
6 1	Fehlersuche	103
		102
6.1	Ubersicht	
6.2	Meldungen auf dem Geratedisplay	104
7 \$	Service	111
7.1	Allgemeine Hinweise und Sicherheitsmaßnahmen	111
7.2	Leaksensor	
7.3	Wechseln der Spritze	
7.4	Wechseln der Pufferschleife	116
7.5	Nadelpaar	118
7	7.5.1 Überprüfen der Nadelpaar-Position	119
7	7.5.2 Waschen der Probennadel	
7	7.5.3 Wechseln der Probennadel	
7	7.5.4 Wechseln des Vorstechers	
7.6	Wechseln der Probenschleife	
7.7	Injektionsventil und Fraktionierventil	
7	7.7.1 Reinigen des Ventils	
7	7.7.2 Auseinanderbauen des Ventils	
7	7.7.3 Zusammenbauen des Ventils	
7.8	Wechseln der Sicherungen	
7.9	Aktualisieren der Autosamplerfirmware	
8	Autosampler-Spezifische Informationen	135
8.1	Funktionsprinzip WPS-3000PL (RS)	
8.2	Autosampler zur Fraktionssammlung	
8	3.2.1 WPS-3000FC	
8	3.2.2 Autosampler mit Micro Fraction Collection (MFC) Option	145
8.3	Biokompatible Autosampler	146
8.4	Injektionsmethoden	150
8	3.4.1 Full-Loop-Injektionen	150
8	3.4.2 Partial-Loop-Injektionen	154
8	3.4.3 Mikroliter Pick-up	157
8	3.4.4 Low Dispersion-Modus	160
9 '	Technische Daten	
10		
10 2	Lubehor, Ersatzteile und Verbrauchsmaterialien	
10.	1 Standardzubehör	
10.	2 Optionales Zubehör.	
1	10.2.1 WPS-3000(T)PL (RS)	171
1	10.2.2 WPS-3000TBPL Nano/Cap	

11 Index	
10.3 Ersatzteile und Verbrauchsmaterialien	
10.2.5 WPS-3000TBFC	
10.2.4 WPS-3000TFC	

# 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> Autosamplers benötigen. Dennoch sollten Sie, bevor Sie mit dem Autosampler arbeiten, die gesamte Anleitung einmal gründlich durchlesen, um sich einen Überblick zu verschaffen.

Alle Beschreibungen innerhalb des Handbuchs beziehen sich auf die folgenden Modelle innerhalb der UltiMate<sup>TM</sup> 3000-Autosamplerserie:

- WPS-3000PL (RS) (einschließlich "T"-Modell mit Probenthermostatisierung)
- WPS-3000TBPL
- WPS-3000TFC
- WPS-3000TBFC

Für die Beschreibungen innerhalb dieses Handbuchs gelten die folgenden Konventionen:

- Für die Beschreibung wird der Ausdruck "der Autosampler" oder "das Gerät" 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-3000PL (RS) gleichermaßen für die Versionen mit Probentemperier-Option (WPS-3000TPL RS).
  - ◆ Viper<sup>™</sup>-Kapillarverbindungen ebenso für nanoViper<sup>™</sup>- und gegebenenfalls andere Viper-Kapillarverbindungen. Teile aus nanoViper bestehen aus PEEK-sheathed Fused Silica, wenn nicht anders angegeben.
- 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 Firmware-Version 4.13 und Chromeleon<sup>™</sup> 6.80 Service Release 13. Wenn Sie den Autosampler unter Chromeleon 7 betreiben möchten, beachten Sie bitte die Hinweise auf Seite 29.

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 diesen Unterlagen 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 Inc. 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.

PEEK ist ein Warenzeichen von Victrex PLC.

PEEKsil ist ein Warenzeichen von SGE Incorporated.

Cheminert ist ein eingetragenes Warenzeichen von Valco Instruments Co., Inc.

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 Sicherheit

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

### 1.2.1 Symbole am Gerät 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 (O)		
<b>1 (1)</b>	Kennzeichnung entsprechend der Richtlinie "Measures for Administration of the Pollution Control of Electronic Information Products" (China-RoHS)		
	Handverletzungsgefahr		
$\triangle$	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.

**I** Hinweis: Bevor Sie mit dem Autosampler beginnen zu arbeiten, lesen Sie diese Anleitung vollständig durch, 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.

#### WICHTIGE SICHERHEITSHINWEISE ZUM WPS-3000TFC UND WPS-3000TBFC AUTOSAMPLER!

Die WPS-3000TFC und WPS-3000TBFC Autosampler wurden AUSSCHLIESSLICH für multidimensionale Trennungen von Biomolekülen mit wässrigen und polaren organischen mobilen Phasen, zum Beispiel Methanol, entwickelt, die üblicherweise für Reversed-Phase (RP) Flüssigchromatographie eingesetzt werden.

#### DAHER GILT: BETREIBEN SIE DIESE BEIDEN AUTOSAMPLER NICHT MIT NICHT-LEITENDEN, UNPOLAREN NORMALPHASEN-ELUENTEN! HIERDURCH KANN EINE FEUER- UND/ODER EXPLOSIONSGEFAHR ENTSTEHEN!

Beachten Sie, dass ein Normalphasen-Kit (Best.-Nr. 6820.0060) NUR für den WPS-3000TFC Autosampler (Best.-Nr. 5824.0020) zur Verfügung steht. Verwenden Sie Normalphasen-Eluenten NUR mit dem WPS-3000TFC Autosampler, nachdem der Autosampler mit dem Normalphasen-Kit umgerüstet wurde. DURCH EINEN AUTOSAMPLER-BETRIEB OHNE DAS ENTSPRECHENDE KIT KANN EINE FEUER- UND/ODER EXPLOSIONSGEFAHR ENTSTEHEN!

Führen Sie KEINE Normalphasen-Anwendungen ohne ordnungsgemäßen Erdungsschutz der Nadel durch. Nicht-leitende, nicht-polare Normalphasen-Eluenten, zum Beispiel Hexan, können zu einer elektrostatischen Aufladung an einem nicht-geerdeten Ventil und einer nicht-geerdeten Nadel führen. DURCH DIE ELEKTROSTATISCHE ENTLADUNG KANN EINE FEUER- UND/ODER EXPLOSIONSGEFAHR ENTSTEHEN. Der WPS-3000TBFC Autosampler (Best.-Nr. 5825.0020) darf NICHT für Normalphasen-Anwendungen eingesetzt werden. Für diesen Autosampler steht kein Umrüstkit zur Verfügung. Verwenden Sie den WPS-3000TBFC Autosampler NIEMALS mit Normalphasen-Eluenten! ANSONSTEN KANN EINE FEUER- UND/ODER EXPLOSIONSGEFAHR ENTSTEHEN. Der WPS-3000TBFC Autosampler darf NUR für Reversed-Phase-Chromatographie eingesetzt werden.

#### • 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>TM</sup> zum Einsatz. Dieses Polymer weist eine ausgezeichnete Beständigkeit gegen die meisten organischen Lösungsmittel auf. Es neigt jedoch dazu aufzuquellen, wenn es mit Trichlormethan (CHCl3), 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.
- 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.
- Ersetzen Sie durchgebrannte Sicherungen immer durch die von Thermo Fisher Scientific autorisierten Original-Ersatzsicherungen.
- Tauschen Sie beschädigte Kommunikationskabel aus.
- Tauschen Sie beschädigte Netzkabel aus. Verwenden Sie nur die für das Gerät bereitgestellten Netzkabel.
- Verwenden Sie ausschließlich die von Thermo Fisher Scientific für das Gerät autorisierten Original-Ersatz- und Zubehörteile.

- 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 Betrieb Ihres HPLC-Systems stets darauf, dass die Minimum-Druckabschaltung gesetzt ist. So vermeiden Sie, dass durch eventuelle Undichtigkeiten Schäden entstehen. Außerdem wird dadurch ein Trockenlaufen der Pumpe verhindert.
- 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.
- Spülen Sie Peroxide bildende Lösungsmittel und Pufferlösungen nach Arbeitsende aus.
- Spülen Sie bei der Umstellung des Lösungsmittels von Puffer auf organische Lösungsmittel das HPLC-System zuvor gründlich mit entionisiertem Wasser oder mit Wasser in HPLC-Qualität.
- Wenn Sie auf ein anderes Laufmittel umstellen, achten Sie auf die Mischbarkeit des neuen Laufmittels mit dem Laufmittel, das im HPLC-System enthalten ist. Sind die Laufmittel nicht mischbar, kann das System beschädigt werden, z.B. durch Ausflockungen.
- Wenn eine Undichtigkeit auftritt, schalten Sie den Autosampler sofort aus und beheben Sie die Ursache für die Undichtigkeit.
- Verwenden Sie nur handelsübliche Lösungsmittel in HPLC-Qualität und Puffer, die kompatibel mit den medienberührten Teilen des Detektors sind.
- Schauen Sie niemals direkt in die LED für die Innenbeleuchtung und verwenden Sie keine lichtbündelnden Instrumente zur Betrachtung des Lichtstrahls. Die hohe Leuchtkraft der Lampe kann schädlich für die Augen sein.
- Um Verletzungen zu vermeiden, greifen Sie bitte während einer laufenden Analyse niemals in den Probenraum.
- Schalten Sie das Ventil des WPS-3000(T)PL RS nicht trocken. Schäden an den Ventilen können die Folge sein.
- Beachten Sie bei längeren Betriebsunterbrechungen (= mehrere Tage) oder wenn Sie den Autosampler zum Versand vorbereiten, die Hinweise zur Außerbetriebnahme des Autosamplers (→ Seite 97).
- Setzen Sie den Autosampler nur entsprechend der bestimmungsgemäßen Nutzung 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 Gerätes

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 für Laborforschungsaufgaben in der HPLC und UHPLC (Ultra-High Performance Liquid Chromatography) entwickelt. 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 kann über das Chromatographie-Management-System Chromeleon gesteuert werden. Als Teil des UltiMate 3000-Systems kann er auch mit anderen Datensystemen betrieben werden, wie

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

Weitere Fragen beantwortet Ihnen gern die Thermo Fisher Scientific-Vertriebsorganisation für Dionex HPLC-Produkte.

Beachten Sie für die Verwendung des Autosamplers Folgendes:

- Die WPS-3000TFC und WPS-3000TBFC Autosampler wurden ausschließlich für multidimensionale Trennungen von Biomolekülen mit wässrigen und polaren organischen mobilen Phasen, zum Beispiel Methanol, entwickelt, die üblicherweise für Reversed-Phase (RP) Flüssigchromatographie eingesetzt werden. Beachten Sie für Normalphasen-Applikationen die Sicherheitshinweise zu den WPS-3000TFC und WPS-3000TBFC Autosamplern auf Seite 5.
- Der Autosampler darf nur mit den original Zubehör- und Ersatzteilen (→ Seite 165) und innerhalb seiner technischen Spezifikationen (→ Seite 132) betrieben werden.
- Verwenden Sie nur handelsübliche Lösungsmittel und Wasser in HPLC-Qualität (besser: in LC-MS-Qualität). Wird Wasser aus Wasseraufbereitungsanlagen verwendet, die nicht ordnungsgemäß gewartet wurden, so kann dies durch polymere Verunreinigungen zu einer starken Beschädigung der Säule und schnelleren Verstopfung der Lösungsmittelfilter führen. Achten Sie darauf, dass Lösungsmittel und Puffer kompatibel mit den medienberührten Teilen des Autosamplers sind.
- Achten Sie auf spezifische Eigenschaften der Lösungsmittel, wie Viskosität, Siedepunkt, UV-Absorption (UV/VIS-Detektor) und Brechungsindex (Brechungsindex-Detektor).

- *Autosampler in Edelstahlausführung* Pufferkonzentrationen: Typischerweise bis zu 1 mol/L. (< 0,1 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.

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.

# 2 Überblick

## 2.1 Kurzbeschreibung

Der Autosampler ist ein hochwertiger Baustein modularer HPLC-Anlagen und Teil des UltiMate 3000-Systems. Er wurde speziell für Anwendungen entwickelt, bei denen hohe Verfügbarkeit und Reproduzierbarkeit auch bei kleinsten Injektionsvolumina wichtig sind. 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 24), die in ein und derselben Autosampler-Sequenz verwendet werden können.
- Die folgenden Injektionsmethoden werden unterstützt: Full-Loop, Partial-Loop und Mikroliter-Pickup. Zur Optimierung des Injektionsprofils bei Full-Loop- und Partial-Loop-Injektionen wird eine Injektionsregelung entsprechend dem "Low Dispersion"-Prinzip unterstützt (→ Seite 150).
- Eine separate Nadel (Vorstecher) durchsticht vor der Injektion die Kappe/das Septum des Probenfläschchens bzw. die Versiegelung der Platte. Damit können Fused-Silica-Kapillaren sicher in den Probenbehälter eingeführt werden. So können selbst kleinste Probenvolumina ohne Probenverlust analysiert werden.
- 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.
- Zur Fehlererkennung stehen im Hinblick auf die Geräteleistung und Zuverlässigkeit diverse Sicherheits- und Überwachungsfunktionen zur Verfügung (→ Seite 22).
- 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 besitzen.
- Der Autosampler steht auch in biokompatibler Ausführung (→ Seite 14) sowie mit integrierter Temperaturregelung zur Kühlung/Erwärmung der Proben (→ Seite 15) zur Verfügung. Darüber hinaus stehen Autosamplerversionen für die Fraktionssammlung zur Verfügung (→ Seite 15).
- Der Probenraum des Autosamplers kann mit Hilfe einer optional verfügbaren transparenten Frontabdeckung (Best.-Nr. 6820.1427) verschlossen werden, z.B. zum Schutz gegen Staub oder andere Partikel.

## 2.2 Funktionsprinzip

Der Autosampler arbeitet nach dem Pulled-Loop Injektionsprinzip und unterstützt die folgenden Injektionsmethoden Full-Loop, Partial-Loop und Mikroliter-Pickup. Zur Optimierung des Injektionsprofils bei Full-Loop- und Partial-Loop-Injektionen wird eine Injektionsregelung entsprechend dem "Low Dispersion"-Prinzip unterstützt.

Eine schematische Darstellung der Funktionsprinzipien sowie detaillierte Informationen zu den Injektionsmethoden finden Sie im Kapitel 8 ( $\rightarrow$  Seite 135).

Folgende Informationen finden Sie	Auf Seite
Funktionsprinzip WPS-3000PL (RS)	135
Funktionsprinzip WPS-3000FC	138
Injektionsmethoden - Full-Loop Injektion - Partial-Loop Injektion - Mikroliter-Pickup - Low Dispersion Modus	150 150 154 157 160

## 2.3 Autosamplerkonfigurationen

## 2.3.1 Übersicht

Der Autosampler ist in den in den Tabellen genannten Konfigurationen verfügbar. Informationen zu den optional erhältlichen Komponenten finden Sie auf Seite 171.

WPS-3000PL RS	BestNr.
<ul> <li>WPS-3000PL RS Nano/Cap (ohne Probenthermostatisierung) mit</li> <li>2,4 μL-Probennadel</li> <li>25 μL-Spritze</li> <li>50 μL-Pufferschleife</li> <li>1 μL-Probenschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet für einen Druck &lt; 90 MPa (12500 psi)</li> </ul>	5826.0010
WPS-3000TPL RS Nano/Cap, wie WPS-3000PL RS Nano/Cap, jedoch mit Probenthermostatisierung	5826.0020
Zur Durchführung von biokompatiblen Anwendungen mit einem WPS-3000(T)PL RS Autosampler ist das PAEK Modification Kit erhältlich. Das Kit enthält alle Komponente für die Umrüstung der nicht-biokompatiblen Fluidik eines WPS-3000(T)PL RS Autosamplers zu einem biokompatiblen Autosampler. Für weitergehende Informationen sowie die Umrüstung wenden Sie sich an den Thermo Fisher Scientific Kundendienst für Dionex HPLC Produkte.	

WPS-3000PL	BestNr.
<ul> <li>WPS-3000PL Nano/Cap (ohne Probenthermostatisierung) mit</li> <li>2,4 μL-Probennadel</li> <li>25 μL-Spritze</li> <li>50 μL-Pufferschleife</li> <li>1 μL-Probenschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet für einen Druck &lt; 34 MPa (5000 psi)</li> </ul>	
WPS-3000TPL Nano/Cap, wie WPS-3000PL Nano/Cap, jedoch mit Probenthermostatisierung	
WPS-3000TBPL Nano/Cap (biokompatibler Autosampler), wie WPS-3000PL Nano/Cap, jedoch mit Probenthermostatisierung	

WPS-3000TBPL Analytical	BestNr.
<ul> <li>WPS-3000TBPL Analytical (biokompatibler Autosampler) mit</li> <li>15 μL-Probennadel</li> <li>100 μL-Spritze</li> <li>500 μL-Pufferschleife</li> <li>50 μL-Probenschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) geeignet für einen Druck &lt; 34 MPa (5000 psi)</li> <li>Probenthermostatisierung</li> </ul>	5823.0020

WPS-3000FC	BestNr.	
<ul> <li>WPS-3000TFC mit <ul> <li>15 μL-Probennadel</li> <li>250 μL-Spritze</li> <li>500 μL-Pufferschleife</li> <li>50 μL Probenschleife</li> <li>Injektionsventil (2 Positionen, 6 Ports) <ul> <li>geeignet f ür einen Druck &lt; 34 MPa (5000 psi)</li> </ul> </li> <li>Fraktionierventil (2 Positionen, 6 Ports) <ul> <li>geeignet f ür einen Druck &lt; 3,4 MPa (500 psi)</li> </ul> </li> <li>Probenthermostatisierung</li> </ul></li></ul>	5824.0020	
WPS-3000TBFC, wie WPS-3000TFC, jedoch biokompatibler Autosampler	5825.0020	
Beide Autosampler-Versionen können auch für Nano/Kap-Anwendungen optimiert werden (→ Seite 143). Der WPS-3000TFC mit Stahlfluidik kann optional mit einem 90 MPa/900-bar Ventil für RSLC-Anwendungen ausgestattet werden.		
Beachten Sie, dass ein Normalphasen-Kit (BestNr. 6820.0060) nur für den WPS-3000TFC Autosampler (BestNr. 5824.0020) zur Verfügung steht. Verwenden Sie Normalphasen (NP)-Eluenten nur mit dem WPS-3000TFC Autosampler, nachdem der Autosampler mit dem Normalphasen-Kit umgerüstet wurde. Beachten Sie für NP- Applikationen die Sicherheitshinweise zu den WPS-3000TFC und WPS-3000TBFC Autosamplern auf Seite 5.		

## 2.3.2 Biokompatible Autosampler

Nahezu alle Beschreibungen der Standardautosampler (Edelstahl) treffen auch auf die biokompatiblen Versionen zu. Trifft eine bestimmte Beschreibung nur auf eine Version zu, so ist dies entsprechend gekennzeichnet.

Bitte beachten Sie auf jedem Fall die im Kapitel 8.3 aufgeführten Unterschiede am Injektionsventil und den daran angeschlossenen Komponenten ( $\rightarrow$  Seite 146).

## 2.3.3 Autosampler mit Probenthermostatisierung

Der Autosampler ist auch mit integrierter Temperaturregelung zur Probenkühlung bzw. Probenerwärmung erhältlich.

Thermoelektrische (Peltier-) Elemente kühlen oder erwärmen die Kühlflüssigkeit und diese wiederum den Drehteller und die darin befindlichen Komponenten. Diese patentierte Technik erlaubt eine präzise Temperierung der Proben. Die untere Temperaturgrenze für die Kühlung beträgt +4 °C. Darüber hinaus können die Proben auf max. +45 °C erwärmt werden. Ausgehend von der Umgebungstemperatur können die Proben um max. 22 °C gekühlt werden.

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

Bitte beachten Sie die folgenden Hinweise:

- Zur Ableitung von eventuell entstehendem Kondenswasser kann an der Unterseite rechts hinten am Gerät ein Ablaufschlauch angeschlossen werden. Der gekühlte Autosampler verfügt über eine aktive Kondenswasserdrainierung (mittels einer Pumpe).
- Achten Sie beim Anschluss des Schlauches 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.
- Bei hoher Luftfeuchtigkeit und starker Kondenswasserbildung wird empfohlen, den Drehteller von Zeit zu Zeit auszuwischen, um Verunreinigungen zum vermeiden. In diesen Fällen kann auch die Verwendung der transparenten Frontabdeckung (Best.-Nr. 6820.1419) hilfreich sein, um die Kondenswasserbildung zu minimieren.

#### 2.3.4 Autosampler zur Fraktionssammlung

Zur Fraktionssammlung stehen folgende Optionen zur Verfügung:

• WPS-3000TFC

Der Autosampler ist mit einem Injektionsventil (2 Positionen, 6 Ports) und einem Fraktionierventil (2 Positionen, 6 Ports) ausgestattet. Eine schematische Darstellung des Funktionsprinzips und Hinweise, die Sie beim Betrieb des Autosamplers beachten sollten, finden Sie im Kapitel 8.2.1 ( $\rightarrow$  Seite 137).

Beachten Sie, dass für Normalphasen-Anwendungen mit dem WPS-3000TFC Autosampler ein Normalphasen-Kit (Best.-Nr. 6820.0060) zur Verfügung steht. Weitere Informationen finden in Kapitel 8.2.1.5 ( $\rightarrow$  Seite 144). Beachten Sie für Normalphasen-Applikationen die Sicherheitshinweise zu den WPS-3000TFC und WPS-3000TBFC Autosamplern auf Seite 5.

 WPS-3000(T)PL Nano/Cap und WPS-3000TBPL Nano/Cap Bei diesen Autosamplern kann das Injektionsventil mit 2 Positionen und 6 Ports durch ein Ventil mit 2 Positionen und 8 Ports für die Fraktionssammlung ersetzt werden (→ Seite 145).

### 2.3.5 Kombinationsmöglichkeiten von Spritzen und Schleifen

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

Autosampler	Spritze	Pufferschleife	Probenschleife
WPS-3000PL RS Nano/Cap, WPS-TPL RS Nano/Cap, WPS-3000(T)PL Nano/Cap und WPS-3000TBPL Nano/Cap	25 μL 100 μL 250 μL <sup>1</sup>	50 μL 500 μL 500 μL	1 μL, 5 μL, 10 μL, 20 μL 5 μL, 10 μL, 20 μL, 50 μL 50 μL, 125 μL
WPS-3000TBPL Analytisch	100 μL 250 μL	500 μL 1000 μL	5 μL, 10 μL, 20 μL, 50 μL 250 μL <sup>2</sup>
WPS-3000TFC, WPS-3000TBFC	25 μL <sup>3</sup> 100 μL 250 μL 250 μL	50 μL 500 μL 500 μL 1000 μL	1 μL, 5 μL, 10 μL, 20 μL 5 μL, 10 μL, 20 μL. 50 μL 50 μL, 125 μL 250 μL

<sup>1</sup> Unter der Best.-Nr. 6820.0031 bzw. 6821.0031 (für biokompatible Version) sind für die 250 μL-Spritze Kits erhältlich, die bereits eine geeignete Puffer- und Probenschleife (125 μL) enthalten.

<sup>2</sup> Aufgrund des Materials der 250μL-Probenschleife kann das Offset bei der Injektorlinearität für diese Kombination bis zu 11,0 μL betragen.

<sup>3</sup> Nano/Kap Umrüstkits sind für den WPS-3000TFC Autosampler (Best.-Nr. 6824.0030) und den WPS-3000TBFC Autosampler (Best.-Nr. 6825.0030) erhältlich. Sie enthalten alle empfohlenen Teile für eine Nano/Cap-Konfiguration.

Informationen zu den Bestellnummern finden Sie in den Kapiteln 10.2 ( $\rightarrow$  Seite 171) und 10.3 ( $\rightarrow$  Seite 180).

## 2.4 Innenansicht



#### 2.4.1 WPS-3000PL RS und WPS-3000TBPL

Nr.	Beschreibung
1	Frontdeckel (nach oben aufgeklappt)
2	Spritzenventil ( $\rightarrow$ Seite 55)
3	Beleuchtung des Probenraums, hier durch Frontdeckel verdeckt ( $\rightarrow$ Seite 27)
4	Schlauch zur Ableitung des Spritzenabfalls Schlauchverbindung Spritzenventil - Waschport (Versorgung Waschport)
5	Spritze
6	Position für optionale Waschflasche (für 125 mL Waschflüssigkeit; → Seite 52)
7	Menüstift
8	Deckel Drehteller Manuellen Zugriff auf die Proben im Drehteller erhalten Sie, indem Sie den Deckel am schwarzen Griff nach links schwenken. Verschließen Sie den Drehteller, indem Sie den Deckel ganz nach rechts schwenken, bis Kontakt zum Nadelarm besteht. Zum Aufziehen der Probe wird der Deckel dann durch die Bewegungen des Nadelarms automatisch geöffnet und geschlossen.
9	Mikrotiterplatte oder Probenhalter
	Detailansicht Drehteller siehe Abb. 2 ( $\rightarrow$ Seite 18).

Nr.	Beschreibung
10	Wasch-Port Der Wasch-Port befindet sich in einer Auffangwanne, in der durch Undichtigkeiten austretende Flüssigkeiten gesammelt werden. Über den Auslass rechts unterhalb des Autosamplers wird die Flüssigkeit aus dem Waschport in den Abfall abgeleitet (→Seite 56)
11	Leaksensor
12	Auffangwanne
13	Nadelpaar, bestehend aus Vorstecher und Probennadel, mit Sensor automatischen Probenfläschchen- bzw. Mikrotiterplatten- und Deep-Well-Plate- Erkennung
14	Injektionsventil mit Probenschleife ( $\rightarrow$ Seite 28)
15	Pufferschleife ( $\rightarrow$ Seite 28)



Abb. 2: Ansicht Drehteller

Nr.	Beschreibung
1	Mikrotiterplatte oder Probenhalter
2	Halter zur Aufnahme von 5 Reagens- bzw. Transportfläschchen à 10 mL. Drei Halter sind im Drehteller installierbar und werden mitgeliefert.



2.4.2 WPS-3000TFC und WPS-3000TBFC

Abb. 3: Innenansicht (hier WPS-3000TFC)

Nr.	Beschreibung
1	Frontdeckel (nach oben aufgeklappt)
2	Spritzenventil ( $\rightarrow$ Seite 55)
3	Beleuchtung des Probenraums, hier durch Frontdeckel verdeckt ( $\rightarrow$ Seite 27)
4	Schlauch zur Ableitung des Spritzenabfalls Schlauchverbindung Spritzenventil - Waschport (Versorgung Waschport)
5	Spritze
6	Position für optionale Waschflasche (für 125 mL Waschflüssigkeit; $\rightarrow$ Seite 52)
7	Menüstift
8	Deckel Drehteller Manuellen Zugriff auf die Proben im Drehteller erhalten Sie, indem Sie den Deckel am schwarzen Griff nach links schwenken. Verschließen Sie den Drehteller, indem Sie den Deckel ganz nach rechts schwenken, bis Kontakt zum Nadelarm besteht. Zum Aufziehen der Probe wird der Deckel dann durch die Bewegungen des Nadelarms automatisch geöffnet und geschlossen.
9	Mikrotiterplatte oder Probenhalter Detailansicht Drehteller siehe Abb. 2 ( $\rightarrow$ Seite 18).

Nr.	Beschreibung
10	Wasch-Port Der Wasch-Port befindet sich in einer Auffangwanne, in der durch Undichtigkeiten austretende Flüssigkeiten gesammelt werden. Über den Auslass rechts unterhalb des Autosamplers wird die Flüssigkeit aus dem Waschport in den Abfall abgeleitet (→Seite 56)
11	Leaksensor
12	Auffangwanne
13	Nadelpaar, bestehend aus Vorstecher und Probennadel, mit Sensor automatischen Probenfläschchen- bzw. Mikrotiterplatten- und Deep-Well-Plate- Erkennung
14	Fraktionierventil
15	Injektionsventil mit Probenschleife ( $\rightarrow$ Seite 28)

# 

## 2.5 Gerätevorderseite

Abb. 4: Gerätevorderseite (hier: WPS-3000TBFC)

Nr.	Bedienelement	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>Funktionen und Menüs, die über Softkeys aufgerufen werden können (→ Seite 74)</li> <li>Fehlermeldung (→ Seite 103)</li> </ul>
2	Standby- Schalter	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 ändern kann, muss der Standby-Schalter circa 1 Sekunde lang gedrückt werden.
3	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; das System ist in Ordnung. Die LED leuchtet rot, wenn ein Fehler erkannt wurde. Autosampler mit Probentemperierung: Die LED leuchtet gelb, wenn die Solltemperatur noch nicht erreicht ist.

## 2.6 Geräterückseite



Abb. 5: Anschlüsse auf der Geräterückseite (hier: WPS-3000TPL)

Nr.	Beschreibung
1	Netzschalter
2	Sicherungsschlitten ( $\rightarrow$ Seite 23)
3	Netzanschluss ( $\rightarrow$ Seite 35)
4	Typenschild
5	USB 1.1-Schnittstelle (Universal Serial Bus Interface): Anschluss des Autosamplers an den Server-PC ( $\rightarrow$ Seite 34)
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 34)
7	Digital I/O zur Kommunikation mit externen Geräten unter Chromeleon ( $\rightarrow$ Seite 35)
8	Nur Autosampler mit Probenthermostatisierung Ventilatoren

### 2.6.1 Netzschalter

Der Netzschalter befindet sich auf der Geräterückseite. Schalten Sie den Autosampler über diesen Schalter ein. Im Normalbetrieb brauchen Sie den Autosampler nicht über den Netzschalter auszuschalten. Verwenden Sie stattdessen die Standby-Taste auf der Gerätevorderseite (→ Seite 21). Drücken Sie die Taste ca. 1 Sekunde lang, damit der Autosampler den Modus ändern kann. Schalten Sie den Autosampler aber über den Netzschalter aus, wenn Sie dazu aufgefordert werden, zum Beispiel, für bestimmte Wartungsarbeiten oder bei längeren Betriebspausen (Stillstand). Beachten Sie dazu auch die Hinweise auf Seite 97.

#### 2.6.2 Sicherungsschlitten

Im Sicherungsschlitten befinden sich zwei Sicherungen. Abhängig vom Autosampler-Modell sind dies zwei Sicherungen mit 2 A, 250 V (Autosampler ohne Probenthermostatisierung) bzw. 4 A, 250 V (Autosampler *mit* Probenthermostatisierung). Informationen zum Sicherungswechsel finden Sie auf Seite 131.

#### 2.6.3 USB-Anschluss

Der Autosampler kann über eine USB-Verbindung (USB = Universal Serial Bus) über das Chromatographie-Management-System **Chromeleon** gesteuert werden. Die Datenübertragung erfolgt dabei digital über das entsprechende USB-Anschlusskabel ( $\rightarrow$  Seite 34). Verwenden Sie zum Anschluss nur die mitgelieferten Anschlusskabel. Nur so wird eine einwandfreie Funktion der Verbindung gewährleistet.

Der interne USB-Hub ( $\rightarrow$  Abb. 5, Nr. 6) erlaubt den Anschluss von drei weiteren Geräten der UltiMate 3000-Reihe an den Autosampler bzw. eines externen USB-Hubs.

**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 Kapitel 3.4.1 ( $\rightarrow$  Seite 34).

## 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. Dazu ist ein entsprechendes Anschlusskabel erforderlich (Best.-Nr. 6000.1004).

Informationen zum Anschluss finden Sie auf Seite 35.

## 2.7 Drehteller, Probenhalter, Probengefäße 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 58.



Abb. 6: Probenhalter (Beispiele)

Folgende Probenhalter stehen für Probenfläschchen zur Verfügung:

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

Wenn Sie Mikrotiterplatten verwenden, müssen Sie diese ebenfalls auf einen geeigneten Probenhalter aufsetzen. Folgende Probenhalter stehen für Mikrotiterplatten zur Verfügung:

Probenhalter für	BestNr.
Standard-Mikrotiterplatten (12 - 24 mm Höhe)	6820.4070, 6820.4086, 6820.4087, 6820.4090 oder 6820.4091
Deep-Well-Plates (34 - 46 mm Höhe)	6820.4079
Deep-Well-Plates (30 - 36 mm Höhe)	6820.4083
Deep-Well-Plates (20 - 32 mm Höhe)	6820.4089
Low-Well-PCR-Plates (8 - 12 mm Höhe)	6820.4070, 6820.4086, 6820.4087, 6820.4090 oder 6820.4091
	Zusätzlich ist ein Adapter (BestNr. 6820.4088) erforderlich.

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 61).

Informationen zu den von Thermo Fisher Scientific empfohlenen Mikrotiterplatten und Probenfläschchen sowie zu den erforderlichen Gefäßhöhen entnehmen Sie bitte der folgenden Tabelle. 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-Vertriebsorganisation für Dionex HPLC-Produkte.

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
1,1-mL-Probenfläschchen (31 - 43 mm)	Konisches Probenfläschchen mit Bördelkappe und geschlitztem Silikon/PTFE Septum für 40-Positionen-Probenhalter (BestNr. 6820.4087)	6000.0077 6000.0076
1,1-mL-Probenfläschchen (31 - 43 mm)	Konisches Probenfläschchen mit Schraubkappe und Silikondichtung für 40-Positionen-Probenhalter (BestNr. 6820.4087)	6000.0078 6000.0057 6000.0058
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
1,8 mL-Probenfläschchen (31 - 43 mm)	Zylindrisches Probenfläschchen mit Bördelkappe und Septum für 40-Positionen-Probenhalter (BestNr. 6820.4070)	6000.0072 6000.0071

Probengefäß (Min Max. Höhe)	Beschreibung	BestNr.
2 mL-Probenfläschchen (31 - 43 mm)	Zylindrisches Probenfläschchen mit Schraubkappe und Silikondichtung für 40-Positionen-Probenhalter (BestNr. 6820.4070)	6000.0060 6000.0057 6000.0058
4-mL-Probenfläschchen (36 - 48 mm)	Zylindrisches Probenfläschchen mit 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
1,5 mL-Eppendorf-Gefäß	Eppendorf-Gefäße ohne Deckel oder mit Deckel, wenn diese als durchstechbar spezifiziert sind; z.B. Eppendorf Standard Reaktionsgefäß 3810; für 40-Positionen Eppendorf-Halter (BestNr. 6820.4094)	Nicht bei Thermo Scientific erhältlich
0,5 mL-Eppendorf-Gefäß	Eppendorf-Gefäße ohne Deckel oder mit Deckel, wenn diese als durchstechbar spezifiziert sind; für 40-Positionen Eppendorf-Halter (BestNr. 6820.4096)	Nicht bei Thermo Scientific erhältlich
Standard-Mikrotiterplatte (12 - 24 mm)	<ul> <li>96 Kavitäten</li> <li>384 Kavitäten</li> <li><i>Hinweis:</i> Setzen Sie diese Platten auf einen der auf</li> <li>Seite 24 genannten Probenhalter auf.</li> <li>Verschlussmatte für Mikrotiterplatte mit</li> <li>96 Kavitäten</li> <li>384 Kavitäten</li> </ul>	6820.4100 6820.4110 6820.4102 6820.4112
Deep-Well-Plates (34 - 46 mm)	96 Kavitäten Verschlussmatte für Mikrotiterplatte mit 96 Kavitäten 384 Kavitäten 24 Kavitäten	6820.4101 6820.4102 Auf Anfrage Auf Anfrage
	<i>Hinweis:</i> Für alle Deep-Well-Plates wird ein spezieller Plattenhalter benötigt (siehe Tabelle Seite 24). Installieren Sie Deep-Well-Plates <i>nicht</i> auf einem der anderen Probenhalter.	
	Für die 24-Deep-Well-Plates ist ein Adapter für den Niederhalter erforderlich (BestNr. 6820.2402).	

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. Wird der Adapter installiert, *muss* der automatische Tray-Test ausgeschaltet werden (→ Seite 95).
- 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 Autosamplerbetriebs 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 dauerhaft aktiviert. Setzen Sie gegebenenfalls das Property **TrayIllumination** in Chromeleon auf **Off**, wenn Ihre Proben lichtempfindlich sind.

## 2.9 Fluidische Anschlüsse

Die fluidischen Anschlüsse im Autosampler sind leicht zugänglich ( $\rightarrow$  Seite 17). Klappen Sie dazu 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- Ersatzkapillaren.
	Bereits benutzte Fittingverbindungen mit Schneidringen aus Stahl sollten nur für dieselbe Kapillarverbindung wieder verwendet werden, um ein erhöhtes Totvolumen zu vermeiden. nanoViper- Kapillaren können mehrmals auch für unterschiedliche Verbindungen benutzt werden.

## 2.10 Leaksensor

Der Autosampler verfügt über einen Leaksensor. Wenn sich Flüssigkeit in der Auffangwanne sammelt, spricht der Leaksensor an und die LED **Status** auf der Gerätevorderseite leuchtet rot. Je nach **Leak sensor mode** erscheint außerdem eine Meldung auf dem Gerätedisplay und es ertönt ein akustisches Signal ( $\rightarrow$  Seite 81).

Wenn der Leaksensor angesprochen hat, beseitigen Sie die Undichtigkeit und trocknen Sie den Leaksensor ( $\rightarrow$  Seite **Fehler! Textmarke nicht definiert.**). Die LED **Status** bleibt rot, bis der Sensor trocken ist. Wählen Sie **Clear** auf der Navigationsleiste, um die Meldung auf dem Gerätedisplay zu löschen ( $\rightarrow$  Seite 78).

## 2.11 Injektionsventil

Das Injektionsventil (2 Positionen, 6 Ports, Cheminert<sup>®</sup>) im Autosampler ist für Anwendungen mit folgendem Druck geeignet:

Injektionsventil	Material	Maximaler Druck	Beschreibung
WPS-3000PL RS	Edelstahl	90 MPa (900 bar)	1/16", 0,25 mm, 50°C
WPS-3000PL, WPS-3000TFC	Edelstahl	35 MPa (350 bar)	1/16", 0,25 mm, 75°C
WPS-3000TBPL, WPS-3000TBFC	PAEK	35 MPa (350 bar)	1/16", 0,25 mm, 50°C

Das Injektionsventil hat zwei Betriebspositionen: Load und Inject. In der Position Load wird die Probe über die Probennadel in die Probenschleife gesaugt. Wenn das Injektionsventil in die Position Inject geschaltet wird, wird die Probe in den Hochdruckflussweg zur Analyse auf die Trennsäule aufgebracht.

Die Komponenten sind wie folgt am Ventil angeschlossen:



Abb. 7: Anschlüsse am Injektionsventil (hier: WPS-3000PL RS)

Port	Anschluss von		Port	Anschluss von
1	Spritze/Pufferschleife		4	Trennsäule
2	Probenschleife		5	Probenschleife
3	Pumpe		6	Probennadel

Wenn Sie mit einem biokompatiblen Autosampler arbeiten, beachten Sie bitte die Hinweise zum Anschluss der Kapillaren am Injektionsventil ( $\rightarrow$  Seite 146).

Der WPS-3000TFC verfügt über *zwei* Schaltventile, Injektions- und Fraktionierventil. Informationen zum Fraktionierventil finden Sie im Kapitel 8.2.1.2 ( $\rightarrow$  Seite 139).
# 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**. Für den WPS-3000TFC ist zusätzlich eine Fraktionierlizenz erforderlich ( $\rightarrow$  Seite 141).

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) oder auf einem Steuerfenster (Control Panel). Direkte Befehle werden mit der Eingabe ausgeführt. 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 Hilfe eines Software-Assistenten erstellen oder manuell, indem Sie ein vorhandenes Programm editieren. Weitere Informationen zum automatischen Betrieb finden Sie auf Seite 71.

# **i** 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.

Bitte beachten Sie auch, dass

- Chromeleon 7 eine andere Terminologie verwendet als Chromeleon 6.80. Informationen hierzu finden Sie in dem Dokument 'Glossary -Chromeleon 7.0' (enthalten im Ordner 'Documents' der Chromeleon 7-Installation).
- in Chromeleon 7 gegebenenfalls noch nicht alle Funktionen von Chromeleon 6.80 unterstützt werden.

# 2.13 Wellness und Predictive Performance

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

- Interne Überwachung aller mechanischen Operationen
- Automatischer Selbsttest beim Einschalten
- Automatischer Überlastungsschutz der Probennadel (Vorstecher)
- Automatische Mikrotiterplatten- und Probenflaschenerkennung ( $\rightarrow$  Seite 95)
- Leaksensor ( $\rightarrow$  Seite 112)
- 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 sowie Service- und (Re)qualifizierungsmaßnahmen zur Verfügung, mit denen ein rechtzeitiger Austausch sichergestellt werden kann (= Predictive Performance;  $\rightarrow$  Seite 91).

# 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 132) 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 stabil sein.
- Vermeiden Sie direkte Sonneneinstrahlung und hohe Luftfeuchtigkeit.
- Achten Sie bei der Aufstellung des Gerätes 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 überprüfen Sie bitte 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, da nur bei sofortiger Reklamation die Transportversicherung für die aufgetretenen Schäden aufkommt.

> 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 Sortimentskasten mit dem Zubehör sowie das Netzkabel. Die beweglichen Teile für den Drehteller (Probenhalter, Mikrotiterplatte) und die Waschflasche sind separat verpackt.

- 2. Heben Sie das Gerät vorsichtig aus dem Karton, und stellen Sie es auf eine stabile Unterlage.
  - Vorsicht: Der Autosampler wiegt mehr als 20 kg. Daher sollten Sie ihn nicht allein anheben. Um ein Herunterfallen des Gerätes zu vermeiden, greifen Sie bitte den Autosampler selbst an der Geräteunterseite, nicht die seitlichen Verpackungsteile. Wenn Sie den Autosampler anheben oder bewegen möchten, greifen Sie bitte immer unter den Boden oder heben Sie das Gerät an den Seiten an.
- 3. Entfernen Sie nun die Verpackungsteile und die Kunststoff-Folie.
- 4. Entfernen Sie die Verpackungsschaumteile, die den Autosampler während des Transports schützen:
  - Zwei Schaumteile arretieren den Frontdeckel. Klappen Sie den Frontdeckel nach oben auf und entnehmen Sie die beiden Schaumteile.
  - Der Nadelarm und Drehteller werden durch ein Verpackungsschaumteil gesichert, das auch als Transportverpackung für die Kapillare zwischen der Pumpe und dem Autosampler dient. Entfernen Sie das Schaumteil und entnehmen die Kapillare.
- 5. Temperieren Sie den Autosampler nach dem Auspacken mindestens vier Stunden, damit sich kondensierter Niederschlag (Betauung), der sich eventuell während des Transports gebildet hat, verflüchtigen kann. Schließen Sie den Autosampler dabei nicht an die Netzstromversorgung an. Bestehen offensichtliche Zweifel an der vollständigen Auflösung des Niederschlags, ist der Autosampler so lange im nicht angeschlossenen Zustand (Netzstecker ausgesteckt) zu akklimatisieren, bis sich der Niederschlag vollständig verflüchtigt hat.

# 3.3 Position des Autosamplers im UltiMate 3000-System

Wird der Autosampler innerhalb eines UltiMate 3000-Systems betrieben, wird empfohlen, die Module des Systems, zum Beispiel wie in Abb. 8 übereinander anzuordnen. Der individuelle Systemaufbau hängt jedoch von der jeweiligen Anwendung ab.



Abb. 8: Beispiel für die Aufstellung eines UltiMate 3000 RSLCnano-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, sollten Sie sich vergewissern, 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 und 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 auf der Geräterückseite mit dem Rechner ( $\rightarrow$  Abb. 5, Seite 22). Wählen Sie eine der folgenden Alternativen:

- Verbinden Sie den Autosampler direkt mit dem USB-Anschluss des Rechners.
- Verbinden Sie den Autosampler mit dem internen USB-Anschluss eines anderen Moduls innerhalb des UltiMate 3000-Systems, das mit dem Rechner verbunden ist.

**Hinweis:** Genauere Informationen dazu, wie die Geräte auf der Rückseite miteinander verbunden werden, finden Sie in weiteren Dokumenten:

- WPS-3000(T)PL RS: im Handbuch zum UltiMate 3000 Nano-Cap-System NCS
- WPS-3000(T)PL, WPS-3000TBPL Nano/Cap: im UltiMate 3000 Proteomics MDLC-Systemhandbuch
- WPS-3000TBPL Analytisch: im Handbuch zur UltiMate 3000 Pumpe
- Verbinden Sie den Autosampler über einen externen USB-Hub mit dem Rechner.

Für den Anschluss stehen die folgenden Kabel zur Verfügung:

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 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 der Geräterückseite ( $\rightarrow$  Abb. 5, Seite 22) das 6-adrige Mini-DIN-Signalkabel (Best.-Nr. 6000.1004).

- 1. Verbinden Sie den 6-poligen Stecker des Kabels mit dem Port **Digital I/O 1 (2, 3** oder **4**).
- Verbinden Sie f
  ür jeden Relaisausgang und digitalen Eingang, den Sie verwenden m
  öchten, die Ader des Kabels mit dem entsprechenden Anschluss des externen Ger
  ätes (→ Dokumentation f
  ür das externe Ger
  ät).
- 3. Nehmen Sie bei der Konfiguration des Autosamplers im Programm **Server Configuration** von Chromeleon (→ Seite 38) folgende Einstellungen vor:
  - Wählen Sie auf der Seite **Relays** den entsprechenden Relaisausgang aus ( $\rightarrow$  Seite 46).
  - ♦ Wählen Sie auf der Seite Inputs den entsprechenden digitalen Eingang aus (→ Seite 46).

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. 9: 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. 10: Digital I/O

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:

Pin	Signalname	Signalpegel	Farbe Ader	Label Ader
1	nicht belegt		pink	Analog High
2	Relay_NC	max. 100VDC/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. 100VDC/0,25A/3W	braun	Digital Output
8	Relay_COM	max. 100VDC/0,25A/3W	weiß	Dig./Out (GND)
			schwarz	Shield

Abb. 11: Anschlussbelegung (6-adrige Mini-DIN Buchse und Kabel)

#### Synchronisation der Injektion mit einem Massenspektrometer

Relay\_4 wird vom Autosampler verwendet, um ein Inject-Response-Signal auszugeben. Dies dient in der Regel dazu, um automatisch vom Autosampler ein Startsignal an ein Massenspektrometer zu senden. Verbinden Sie in diesem Fall das Massenspektrometer mit dem Ausgang Digital I/O 4. Es ist nicht notwendig, das Relay\_4 im Programm Server Configuration zu aktivieren. Die Dauer des Signals kann im Dialogfenster **Commands** in Chromeleon festgelegt werden:

- 1. Drücken Sie F8 oder klicken Sie Command im Menü Control.
- 2. Klicken Sie auf das Pluszeichen neben Sampler.
- 3. Wählen Sie **Relay4Enabled**.
- 4. Wählen Sie eine Dauer zwischen 1 s (**1sec\_InjectOut**) und 10 s (**10sec\_InjectOut**) aus der Liste aus.
- 5. Klicken Sie auf Execute und schließen Sie das Dialogfenster.

Wenn Sie das Massenspektrometer *nicht* sofort nach der Injektion starten möchten, können Sie ein anderes Relais konfigurieren und das Startsignal manuell ausgeben. Das Relais muss im Programm Server Configuration aktiviert und die entsprechenden Relay On/Off-Kommandos in das Programm eingefügt werden.

Hinweis: Wenn Sie Relay\_4 nicht für die Synchronisation der Injektion verwenden möchten, können Sie das Relay\_4 für die benutzerdefinierte Steuerung unter Chromeleon aktivieren, indem Sie in Chromeleon das Property Relay4Enabled auf Yes setzen.

# 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 Hilfe zu Chromeleon.

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

#### 3.5.1 Laden des USB-Treibers für den Autosampler

- 1. Schalten Sie, sofern noch nicht geschehen, den Chromeleon-Rechner ein.
- 2. Melden Sie sich unter Windows<sup>®</sup> 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.
- 3. Starten 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.
- 5. Schließen Sie das Server Monitor-Fenster mit Close. Das Server Monitor-Symbol serscheint in der Taskleiste.

**i** Hinweis: Über den Knopf 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.

Wenn Windows den Autosampler nicht automatisch erkennt und stattdessen einen Installationsassistenten startet, 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 und ziehen Sie das USB-Kabel vom Rechner ab.
- c) Installieren Sie Chromeleon.
- d) Verbinden Sie das USB-Kabel wieder mit dem Rechner und schalten Sie den Autosampler ein. Windows erkennt jetzt den Autosampler und installiert die USB-Software f
  ür den Autosampler automatisch.

#### 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 und ziehen Sie das USB-Kabel vom Rechner ab.
- c) Installieren Sie Chromeleon.
- d) Verbinden Sie das USB-Kabel wieder mit dem Rechner und 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 38), können Sie den Autosampler in Chromeleon installieren und konfigurieren:

- 1. Starten Sie, falls erforderlich, den Chromeleon Server Monitor ( $\rightarrow$  Seite 38).
- 2. 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 entsprechende Meldung mit OK.
- Überprüfen 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 41).
- 9. Klicken Sie auf **Fertigstellen**, 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 47).

**Hinweis:** Änderungen, die Sie im Dialogfenster **Commands**, in einem Programm (PGM) oder in 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		X
Device Name:	Sampler	
Demo Mode	Firmware Download	
Module Address:		
		Browse
1		

Abb. 12: Seite General

#### • Device Name

Name, unter dem der Autosampler in der Installationsumgebung und im Chromeleon Client geführt wird. Wenn Sie den Autosampler über vorhandene Steuerfenster steuern möchten, sollten Sie den vorgegebenen Namen nicht ändern. 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 gewünschten 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 bitte der Beschreibung im Kapitel 7.9 ( $\rightarrow$  Seite 132).

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 das UltiMate 3000-System eine Niederdruckgradientenpumpe aus der UltiMate 3000-Pumpenserie enthält, sollten Sie hier außerdem den Autosampler mit der Pumpe verbinden. Thermo Fisher Scientific empfiehlt, diese Einstellung grundsätzlich vorzunehmen, da sie erlaubt, den Injektionsbefehl des Autosamplers mit den Kolbenhüben einer Niederdruckgradientenpumpe 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	x
_ Well Plate S	Sampler Segment Types	
Red	40 Standard Vials	
Green	40 Standard Vials	
Blue	40 Standard Vials	
Waming: E	Ensure that these settings match the installed items!	
- Pump Link - Flow throug	igh Sampler is delivered by pump (s):	
<none></none>	<b>_</b>	
<none> UM3PUM UM3PUM UM3PUM UM3PUM</none>	IP_L_STRK (DGP-3600SD Pump)           IP_R_STRK (DGP-3600SD Pump)           IP_STROKE (LPG-3400SD Pump)           S_STRK 1 (WPS-3000(RS) Autosampler #4)	
UM3WPS UM3WPS	S_STRK_2 (WPS-3000(RS) Autosampler #4)	

Abb. 13: 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. Überprüfen und ändern Sie diese Einstellungen gegebenenfalls. Wenn Sie zu einem späteren Zeitpunkt andere Probengefäße im Autosampler installieren, müssen Sie die Einstellung in Chromeleon entsprechend anpassen ( $\rightarrow$  Seite 87).

• Pump Link

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

• Wenn das UltiMate 3000-System eine Pumpe des Typs DGP-3600 oder ein Nano/Cap System NCS-3500 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, LPG-3400 oder einem NCS-3500 verbunden haben, ist die Funktion zur Synchronisierung des Injektionsbefehls mit den Kolbenhüben der Pumpe aktiviert. 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 86).

#### Seite Options

Auf dieser Seite wird die Konfiguration des Autosamplers angezeigt. Prüfen und ändern Sie gegebenenfalls die angezeigte Konfiguration oder nehmen Sie zusätzliche Einstellungen vor:

Options				×
Needle Siz © 2.4 µ © 30 µl	Needle Size		Loop Size	С 100 µ С 125 µ С 250 µ С 500 µ
Syringe Si © 25 µl © 500 j	С 100 µl J С 1000 µl	С 250 µl С 2500 µl	С 50 µl С Othersize (µl):	1
<ul> <li>Thermostatting option installed</li> <li>WPS-3000TBPL Analytical</li> <li>Micro Fraction Collection option installed</li> <li>WPS-3000TFC/WPS-3000TBFC</li> </ul>		lled I ption installed )TBFC	− Buffer Tubing Size −	С 1000 µ С 2000 µ 50
Bridge Volume (µl): 0 ✓ Automatically Wash after Abort errors			Waming: Ensure that the setti page match the inst	ngs on this alled items!

Abb. 14: Seite Options

- Needle Size, Syringe Size, Loop Size und Buffer Tubing Size Zeigen an, welche Nadel-, Spritzen-, Schleifen- und Pufferschleifengrößen installiert sind.
- Thermostatting option installed Vergewissern Sie sich, dass das Kontrollkästchen aktiviert ist, wenn es sich bei Ihrem Autosampler um ein Gerät mit Probenthermostatisierung handelt.

#### • WPS-3000TBPL Analytical

Vergewissern Sie sich, dass das Kontrollkästchen aktiviert ist, wenn Sie mit einem WPS-3000TBPL Analytisch (biokompatibler Autosampler für analytische HPLC-Anwendungen) arbeiten.

Ist dieses Kästchen *nicht* aktiviert, stehen gegebenenfalls nicht alle Funktionen für analytische Systeme zur Verfügung und es kann zu Problemen bei der Operational Qualification kommen.

#### • Micro Fraction Collection option installed

Aktivieren Sie das Kontrollkästchen, wenn der Autosampler über die Option Micro Fraction Collection verfügt, also das 6-Port Injektionsventil gegen ein 8-Port-Ventil zum Sammeln von Fraktionen ausgetauscht wurde. Das Kontrollkästchen muss *nicht* ausgewählt werden, wenn ein WPS-3000(T)FC installiert ist.

#### • WPS-3000FC/WPS-3000TFC

Vergewissern Sie sich, dass dieses Kontrollkästchen aktiviert ist, wenn Sie mit einem WPS-3000TFC arbeiten. Das Kontrollkästchen muss *nicht* ausgewählt werden, wenn eine Micro Fraction Collection-Option installiert ist.

#### BridgeVolume (µL)

Wenn die beiden Ventile im WPS-3000TFC für die Fraktionssammlung über eine Kapillare miteinander verbunden sind, *muss* das Volumen der Kapillare im Feld **BridgeVolume** angegeben sein. Vergewissern Sie sich, dass der angegebene Wert mit dem Wert auf dem Label der Kapillare übereinstimmt (6,2  $\mu$ L für analytische bzw. 1  $\mu$ L für Nano-Anwendungen).

Ist das **BridgeVolume** nicht korrekt angegeben, kann es zu Fehlern bei der Injektionssteuerung kommen.

**i** Hinweis: Denken Sie daher daran, den Wert im Feld BridgeVolume entsprechend anzupassen, wenn Sie zwischen den beiden Ventilen eine Kapillare mit einem anderen Volumen installieren oder wenn Sie den Autosampler für eine Applikation einsetzen, bei der die beiden Ventile *nicht* miteinander verbunden sind (in diesem Fall ist BrigdeVolume auf 0 zu setzen).

#### • 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 **Segment / 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 entsprechende Meldung im Chromeleon Audit Trail.

Überprüfen 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 41).

#### **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.

**Hinweis:** Relay 4 wird vom Autosampler verwendet, um ein Inject-Response-Signal auszugeben. Dies dient in der Regel dazu, um automatisch vom Autosampler ein Startsignal an ein Massenspektrometer zu senden  $(\rightarrow$  Seite 37). Wählen Sie daher für die Steuerung unter Chromeleon ein anderes Relais. Wenn Sie Relay 4 verwenden möchten, können Sie Relay 4 aktivieren, indem Sie in Chromeleon das Property Relay4Enabled auf Yes setzen.

#### **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 40).
- 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 41).

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 entsprechend 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** im Menü **File** und schließen Sie das Serverkonfigurationsprogramm.

# 3.6 Einrichten des Autosamplers in DCMSLink

Wenn Sie den Autosampler in DCMSLink einrichten möchten, finden Sie die entsprechenden Informationen im *DCMSLink Installation Guide*, der auf der DCMSLink-DVD im Verzeichnis *Additional Documents*\*DCMSLink User Documents* zur Verfügung steht.

- 1. Installieren und konfigurieren Sie die DCMSLink-Software ( $\rightarrow$  DCMSLink Installation Guide).
- 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 den Schritten in Kapitel  $3.5.2 (\rightarrow$ Seite 40).
- 4. Konfigurieren Sie den Autosampler. Die Vorgehensweise entspricht der Beschreibung in Kapitel 3.5.3 (→ Seite 41).

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

# 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.

Um Schäden am Ventil zu vermeiden, achten Sie beim Betrieb des 900 bar-Ventils (WPS-3000(T)PL RS) darauf, das Ventil nicht trocken zu schalten.

Nachdem Sie den Autosampler wie in den Kapiteln 3.1 bis 3.5 ( $\rightarrow$  Seite 31 und folgende) beschrieben ausgepackt, aufgestellt und angeschlossen haben, bereiten Sie den Autosampler wie folgt für den Betrieb vor:

1. Schließen Sie den Autosampler, je nach Anwendung, fluidisch an die Pumpe sowie an den Flow Manager bzw. den Säulenthermostat oder die Trennsäule an.

Genauere Informationen zu den fluidischen Verbindungen in Nano- und Kap-HPLC-Systemen, finden Sie in weiteren Dokumenten:

- *WPS-3000(T)PL RS:* im Handbuch Standard Applications für UltiMate 3000 RSLCnano-Systeme.
- *WPS-3000(T)PL, WPS-3000TBPL Nano/Cap:* im UltiMate 3000 Proteomics MDLC-Systemhandbuch.
- Alle Autosampler
   Beachten Sie bitte die Hinweise zum Anschluss der Kapillaren am Injektionsventil im Kapitel 4.2 (→ Seite 51).
- *Biokompatible Autosampler* Beachten Sie bitte die Hinweise zum Anschluss der Kapillaren am Injektionsventil im Kapitel 8.3 (→ Seite 146).
- 2. Überprüfen Sie die Anschlüsse am Injektionsventil (→ Seite 28) bzw. am Injektionsund Fraktionierventil (→ Seite 137).
  - ▲ Vorsicht: Es ist wichtig, dass der Inhalt der Probenschleife im Back-Flush-Modus (→ Seite 135) auf die Säule injiziert wird. Achten Sie daher darauf, dass die Anschlüsse der Säule und der Pumpe am Injektionsventil nicht vertauscht werden.

- 3. Installieren und befüllen Sie die Flasche für die Waschflüssigkeit ( $\rightarrow$  Seite 52).
- 4. Überprüfen Sie die Anschlüsse am Spritzenventil ( $\rightarrow$  Seite 55).
- 5. Schließen Sie die Systemdrainage an ( $\rightarrow$  Seite 56).
- 6. Wenn Sie den Autosampler mit Chromeleon betreiben möchten Installieren Sie den Autosampler in Chromeleon ( $\rightarrow$  Seite 38).

Autosampler für die Fraktionssammlung Installieren Sie zusätzlich in Chromeleon den Treiber Fraction Collection ( $\rightarrow$  Seite 141)

- 7. Schalten Sie den Autosampler ein ( $\rightarrow$  Seite 63).
- 8. Spülen Sie die Spritze ( $\rightarrow$  Seite 56).
- 9. Installieren und bestücken Sie die Probenhalter oder Mikrotiterplatten ( $\rightarrow$  Seite 58).
- Bestücken Sie, falls erforderlich, die 5-Positionen Halter f
  ür die 10 mL-Fl
  äschchen
  (→ Seite 61).
- Vergewissern Sie sich, dass das Nadelpaar korrekt in den Wasch-Port f\u00e4hrt (→ Seite 119).
- Nur Autosampler mit Probenthermostatisierung Schalten Sie die Probenthermostatisierung ein und legen Sie gegebenenfalls die Temperatur fest (→ Seite 86).
- 13. Überprüfen und ändern Sie gegebenenfalls die Einstellungen für den Leakerkennung (→ Seite 89).
- Passen Sie, falls erforderlich, die Helligkeit und den Kontrast der Displayanzeige an Ihre Anforderungen an (→ Seite 89).
- 15. Ehe Sie mit der Probenanalyse beginnen, sollten Sie das gesamte System äquilibrieren (→ Seite 62).

# 4.2 Hinweise zum Anschluss von Kapillaren

Beachten Sie beim Anschluss von Kapillaren an den Autosampler bitte 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 daher nur an der dafür vorgesehenen Position.
- Verwenden Sie nur die mitgelieferten, fertig konfektionierten Anschlusskapillaren bzw. Original-Ersatzkapillaren.
- Verwenden Sie möglichst Viper-Kapillarverbindungen. Beachten Sie dabei 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 Hilfe der schwarzen Rändelschraube und *nur* per Hand fest (verwenden Sie *kein* Werkzeug). Die Rändelschraube kann jederzeit leicht von der Kapillare entfernt und wieder aufgesetzt werden.

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.

Kapillaren mit Viper-Fittingen können für unterschiedliche Verbindungen wieder verwendet werden.

• 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, sollten 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.

Bereits benutzte Fittingverbindungen sollten nur für dieselbe Kapillarverbindung wieder verwendet werden, um ein erhöhtes Totvolumen oder Beschädigungen und Undichtigkeiten zu vermeiden.

Lesen Sie zum Anschluss von Fittingverbindungen im biokompatiblen Autosampler auch Kapitel 8.3.

# 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 Racks beinhaltet, können Sie die Waschflüssigkeit online über den Degaser der Pumpe oder des Solvent Racks entgasen.

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

Die Waschflüssigkeit kann wie folgt an den Degaser angeschlossen werden:

- 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 (→ Seite 53)
   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 Autosamplerzubehörs.

**Hinweis:** Um das Entlüften zu beschleunigen kann es hilfreich sein, statt der 25 μLeine 100 μL-Spritze zu verwenden. Dadurch kann sich jedoch die Präzision und Genauigkeit der Injektion verschlechtern.

#### Direkter Anschluss an den Degaser

- 1. Befüllen Sie eine Vorratsflasche mit einer geeigneten Waschflüssigkeit.
- Bereiten Sie zunächst den Ansaugschlauch für die Waschflüssigkeit vor und platzieren 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. 17, Seite 55).
- 5. Kürzen Sie das freie Schlauchende mit Hilfe eines Kapillarschneiders entsprechend Ihren Anforderungen. Achten Sie auf einen rechtwinkligen Schnitt.
- 6. Verbinden Sie das freie Schlauchende mit Hilfe der Rändelschraube und des Fittings aus dem Online Degas Wash Kit mit dem Degaserauslass.
- 7. Entlüften Sie die Spritze, bis keine Luftblasen mehr sichtbar sind ( $\rightarrow$  Seite 56).

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

- 1. Befüllen 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. 17, Seite 55).
- 5. Verbinden Sie das andere Ende dieses Schlauchs mit dem T-Stück aus dem Online Degas Wash Kit.
- 6. Lösen Sie ggf. die Verbindung zwischen dem Degaserauslass und der Pumpe, je nach Pumpentyp am Sternventil oder am Pumpeneinlass, und schließen Sie den vom Degaserauslass kommenden Schlauch 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. 15: Beispiel für den Anschluss der Waschflüssigkeit über die Pumpe

8. Entlüften Sie die Spritze, bis keine Luftblasen mehr sichtbar sind ( $\rightarrow$  Seite 56).

Anschluss des Schlauchs vom Spritzenventil

#### 4.3.2 Anschluss der Waschflasche

Der Autosampler wird mit einer 125 mL-Waschflasche ausgeliefert. Schließen Sie die Waschflasche folgendermaßen an:

- 1. Befüllen 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.
  - **I** Hinweis: Im Hinblick auf bestmögliche Ergebnisse bei isokratischen Trennungen empfehlen wir, als Waschlösung die mobile Phase (ohne Salze oder Modifier) zu verwenden. Für Gradienten empfehlen wir, eine Waschlösung zu verwenden, die dem Startgradienten (ohne Salz oder Modifier) entspricht. Reines Methanol wird als Waschlösung nicht empfohlen.
- 2. Prüfen Sie, ob der Schlauch für die Waschflüssigkeit durch die Öffnung im Deckel ausreichend tief (bis zum Boden) in die Waschflasche geführt ist. (Halten Sie die Flasche ggf. gegen das Licht.)
- Schließen Sie das andere Ende des Schlauchs am Spritzenventil an (→ Abb. 17, Seite 55).
  - **Hinweis:** Um zu verhindern, dass Luft in das System gelangt, sollte der Schlauch beim Anschließen an das Spritzenventil komplett mit der Waschflüssigkeit befüllt sein. Dies können Sie erreichen, indem Sie beispielsweise die Waschflasche zusammendrücken und dabei das Belüftungsloch im Flaschendeckel mit dem Finger zuhalten, um Druck in der Flasche aufbauen zu können.
- 4. Hängen Sie die Halteklammer mit der Waschflasche in den Halter ein. Die Halteklammer muss im Halter einrasten.



ut Halteklammer (Rückansicht) Halter Waschflasche Abb. 16: Einsetzen der Waschflasche

5. Entlüften Sie Spritze, bis keine Luftblasen mehr sichtbar sind ( $\rightarrow$  Seite 56).

Wenn Sie die Waschflasche neu befüllen oder die Waschflüssigkeit austauschen möchten, greifen Sie bitte die Flasche einschließlich der Halteklammer und schieben Sie beides senkrecht nach oben. Die Halteklammer rastet aus. Ziehen Sie die Flasche einschließlich der Halteklammer heraus.

**I** Hinweis: Wenn Sie für Ihre Applikation mehr Waschflüssigkeit benötigen, können Sie alternativ einen längeren Schlauch (mit Flansch für das Ventilfitting) anschließen und eine größere Waschflasche neben den Autosampler stellen. Um diesen längeren Schlauch vollständig zu füllen, müssen Sie den Spülvorgang unter Umständen mehrmals wiederholen.

# 4.4 Anschlüsse am Spritzenventil

Überprüfen Sie am Spritzenventil, ob die Pufferschleife und die Spritze sowie die Schläuche für den Spritzenabfall korrekt angeschlossen sind. Ziehen Sie die Verbindungen gegebenenfalls nach.



Abb. 17: Anschlüsse am Spritzenventil

Wenn Sie die Verbindung der Spritze oder Waschflüssigkeit nachgezogen haben, sollten Sie die Spritze spülen ( $\rightarrow$  Seite 56). Dadurch werden Luftblasen aus der Spritze, dem Spritzenventil und den Ansaugschläuchen für die Reinigungslösung entfernt.

Wenn Sie die Verbindung der Pufferschleife nachgezogen haben, sollten Sie einen Waschzyklus für die Pufferschleife durchführen, um Luftblasen aus der Pufferschleife zu entfernen. Wählen Sie dazu am Gerätedisplay die Funktionstaste **Wash** ( $\rightarrow$  Seite 79), wählen Sie **Wash** im Menü **Control** ( $\rightarrow$  Seite 79) oder führen Sie den Befehl **Wash** in Chromeleon aus.

# 4.5 Anschließen der Systemdrainage

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 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

- Öffnen Sie in Chromeleon das Dialogfenster Commands f
  ür den Autosampler (→ Seite 68).
- 2. Geben Sie unter **PrimeSyringeNumber** 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** ( $\rightarrow$  Seite 74).
- 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).

#### Hinweis zum Troubleshooting

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

- 1. Bauen Sie die Spritze aus ( $\rightarrow$  Seite 113).
- 2. Befüllen 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 wieder ein ( $\rightarrow$  Seite 113).
- 8. Führen Sie einen Waschzyklus durch. Führen Sie dazu den Befehl **Wash** in Chromeleon aus oder wählen Sie am Gerätedisplay die Funktionstaste **Wash** oder wählen Sie **Wash** im Menü **Control**.

## 4.7 Einsetzen und Bestücken eines Probenhalters

Der Autosampler ist so konzipiert, dass Sie die Probenhalter oder Mikrotiterplatten bequem einsetzen können. Eine Tabelle mit den verfügbaren Probenhaltern finden Sie auf Seite 24.



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. Dazu stehen Ihnen folgende Möglichkeiten zur Verfügung:
  - Drehen Sie den Teller per Hand, bis das gewünschte Segment vorn ist.
  - Wählen Sie die Funktionstaste Rotate, bis das gewünschte 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**).
- 3. Wenn Sie eine Mikrotiterplatte verwenden möchten

Setzen Sie die Mikrotiterplatte auf dem Probenhalter an der Position A1 an und drücken Sie sie fest auf den Probenhalter, bis sie einrastet ( $\rightarrow$  Abb. 19).



Abb. 19: Aufsetzen einer Mikrotiterplatte auf einen Probenhalter

4. Setzen Sie den Probenhalter bzw. den Probenhalter mit Mikrotiterplatte wie unten gezeigt in den Drehteller ein. Der Probenhalter muss dabei auf den Positionierstiften einrasten.





Positionierstifte im Drehteller



*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, z.B. zur Aufnahme der Reagensflüssigkeiten, installiert. Wenn Sie *ausnahmsweise* ohne diese Halter arbeiten möchten, achten Sie bitte 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 ganz nach rechts, bis Kontakt zum Nadelarm besteht.

- 7. Wenn der Autosampler über Chromeleon gesteuert wird
  - a) Überprüfen und ändern Sie in Chromeleon gegebenenfalls die Einstellungen für die Probengefäße (→ Seite 87). Die Probennadel kann beschädigt werden, wenn die Einstellungen in Chromeleon nicht mit den tatsächlichen installierten Probengefäßen übereinstimmen.
  - b) Passen Sie gegebenenfalls den Wert f
    ür die Nadelh
    öhe 
    über den Parameter SampleHeight an (→ 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, falls erforderlich, mit den geeigneten 10 mL-Probenfläschchen ( $\rightarrow$  Seite 24).

- Warnung: Um Verletzungen zu vermeiden, greifen Sie während einer laufenden Analyse niemals in den Probenraum.
   Hinweis: Der 5-Positionen-Halter kann alternativ mit Hilfe von Adaptern (Best.-Nr. 6820.4092) mit 2 mL-Probenfläschchen bestückt werden. Setzen Sie bei Verwendung des Adapters den Wert von SamleHeightOffset\_T auf 9mm, um eine Beschädigung der Nadel zu verhindern.
- 1. Schwenken Sie den Deckel des Drehtellers ganz nach links.
- 2. Holen Sie das Segment (rot, grün oder blau) nach vorn, das Sie bestücken möchten:
  - Drehen Sie den Teller per Hand, bis das gewünschte Segment vorn ist.
  - Wählen Sie Funktionstaste Rotate, bis das gewünschte 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 gewünschten 10 mL-Probenfläschchen ( $\rightarrow$  Seite 24).
- 4. Schwenken Sie den Deckel des Drehtellers wieder nach rechts, bis Kontakt zum Nadelarm besteht.
- Wenn der Autosampler über Chromeleon gesteuert wird
   Passen Sie im Dialogfenster Commands (→ Seite 68) gegebenenfalls den Wert für die Nadelhöhe über den Parameter TransLiquidHeight an (→ Seite 93).

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

# 4.9 Äquilibrieren des Systems

Ehe Sie den Autosampler zur Probenanalyse einsetzen, sollten Sie das gesamte UltiMate 3000-System äquilibrieren:

- 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 (z.B. den Säulenofen) auf die für die Anwendung erforderliche Temperatur.
- 3. Stellen Sie die Wellenlänge ein und schalten Sie die Lampen ein. (Die Wellenlängenkalibrierung erfolgt automatisch.)
- 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** (→ Seite 68) oder über die Steuerfenster (→ Seite 69) aus.
- Automatisieren Sie die Äquilibrierung, indem Sie ein Äquilibrierprogramm erstellen und ablaufen lassen (→ Seite 71).
- **Hinweis:** Für den WPS-3000TBPL Analytisch steht Ihnen zusätzlich der **SmartStartup**-Assistent zur Verfügung. 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 gewünschte Temperatur gebracht. Zusätzlich kann SmartStartup die Pumpe automatisch entlüften, die Trennsäule spülen und das HPLC-System äquilibrieren. Wichtige Modulparameter werden automatisch kontrolliert. Weitere Informationen finden Sie in der Hilfe zu Chromeleon.

#### 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 76). 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, zum Beispiel während der Inbetriebnahme oder für Diagnose- oder Wartungsarbeiten, direkt am Autosampler durchgeführt werden können. Dies sind zum Beispiel:

- Ausführen bestimmter Funktionen (Bewegen des Drehtellers)
- Ausführen von Befehlen (Waschen, Anfahren von Servicepositionen)
- Einstellen von Parametern (Leaksensormodus, Solltemperatur für die Probenthermostatisierung)
- Abfragen von Informationen zu Diagnosezwecken
- Abfragen und Ändern der Gerätekonfiguration

Informationen hierzu finden Sie im Kapitel 5.5 ( $\rightarrow$  Seite 74)

### 5.1 Einschalten des Autosamplers

Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite ein:

- Auf dem Gerätedisplay erscheinen kurzzeitig allgemeine Informationen zum Autosampler: Gerätetyp, Seriennummer, Bootloader- und Firmware-Version.
- Der Autosampler führt einen Selbsttest durch. Dabei werden alle wesentlichen Baugruppen auf korrekte Funktion überprüft. 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 entsprechende 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 ( $\rightarrow$  Seite 21). 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 folgende Statusanzeige auf dem Gerätedisplay:

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

Abb. 22: 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 89).
# 5.3 Probenposition

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



Abb. 23: Probenort (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.

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

# Probenposition bei Probenfläschchen, Deep Well-Plates und Mikrotiterplatten

Bei der Angabe der Probenposition steht der Buchstabe für das Segment, in dem sich das Probengefäß befindet, also R, G oder B, an erster Stelle. 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  $(\rightarrow Abb. 24)$ .



Abb. 24: 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 Reagens- und Transportfläschchen steht der Buchstabe für das Segment, in dem sich das Fläschchen befindet, an erster Stelle (R, G oder B), gefolgt von einer Zahl von 1 bis 5.



Abb. 25: 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.
  - **Hinweis:** *Bevor* Sie den Autosampler mit dem Chromeleon-Rechner verbinden und den Autosampler einschalten, sollten Sie sich vergewissern, dass die Chromeleon-Software auf dem Rechner installiert und der Lizenzcode eingegeben ist. Nur dann wird der USB-Treiber für den Autosampler automatisch geladen und das Windows-Betriebssystem kann den Autosampler erkennen, wenn dieser eingeschaltet wird.
- Der Autosampler ist, wie im Kapitel 3.5 beschrieben, in Chromeleon eingerichtet (→ Seite 38).

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** (→ Seite 68) oder auf einem Steuerfenster (Control Panel) (→ Seite 69)
- Automatisch über ein Steuerprogramm (PGM) ( $\rightarrow$  Seite 71)

## 5.4.1 Verbinden mit Chromeleon

- 1. Starten Sie gegebenenfalls den Chromeleon Server Monitor ( $\rightarrow$  Seite 38).
- 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 verschiedene Funktionen zur Lebensdauerprognose von Verschleißteilen zur Verfügung (→ Seite 91).
- 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

- 1. Öffnen Sie ein (beliebiges) Steuerfenster. Die Steuerfenster sind im Chromeleon-Browser in Verzeichnis **Dionex Templates/Panels** abgelegt und können mit einem Doppelklick geöffnet werden.
- Verbinden Sie das Steuerfenster mit der Zeitbasis, auf welcher der Autosampler installiert ist. Wählen Sie dazu 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.

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

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

Welche Parameter und Befehle angezeigt werden, hängt ab von

- ◆ der Chromeleon-Version
- den in der Autosamplerkonfiguration festgelegten Optionen ( $\rightarrow$  Seite 41).
- 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 gewünschten Filter im Menü aus.



Abb. 26: Dialogfenster Commands

6. Vergewissern Sie sich, dass der Autosampler mit Chromeleon verbunden ist. Verbinden Sie den Autosampler gegebenenfalls über den Befehl **Connect**.

Eine Liste aller für den Autosampler verfügbaren Kommandos und Parameter finden Sie in der Hilfe zu Chromeleon. 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. 27, 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 Hilfe zu Chromeleon.

 Klicken Sie auf dem Panel Tabset f
ür Ihre Zeitbasis die Registerkarte Sampler an (→ Abb. 27, Seite 70). 3. Vergewissern Sie sich, dass der Autosampler mit Chromeleon verbunden (connected) ist (die LED der Schaltfläche grün). Verbinden Sie den Autosampler gegebenenfalls über die Schaltfläche **Connect**.



Abb. 27: 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 Hilfe eines 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 die HPLC-Anlage automatisiert herunterzufahren ( $\rightarrow$  Seite 99), oder Programme, um sicherzustellen, dass die Anlage nach einem Stromausfall wie gewünscht weiterarbeitet. Einzelheiten zu diesen Programmen finden Sie in der Hilfe zu Chromeleon.

#### 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. Nehmen Sie auf den einzelnen Seiten des Assistenten die gewünschten Einstellungen vor oder übernehmen Sie die vorgegebenen Werte. Informationen zu den einzelnen Seiten erhalten Sie über die Schaltfläche Hilfe.
- 3. Nach Abschluss des Programm-Assistenten erstellt Chromeleon automatisch das entsprechende Programm.
- 4. Starten Sie das Programm wie unten beschrieben ( $\rightarrow$  Seite 73).

#### Manuelles Erstellen eines Programms

1. Öffnen Sie ein vorhandenes Programm.

Öffnen Sie das gewünschte 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 gewünschte Programm aus.



Abb. 28: Chromeleon Programm (hier: Programmansicht im Commands View)

2. Nehmen Sie in dem Programm die gewünschten Änderungen vor. Die Geräteansichten (Device Views) bieten die einfachste Möglichkeit, ein Programm zu ändern. Klicken Sie ein Gerät an und nehmen Sie die gewünschten 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 Hilfe zu Chromeleon.

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 (Peakidentifizierung, Flächenund Stoffmengenbestimmung) enthalten.
- 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 Hilfe zu Chromeleon.

#### Andere Programme

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

# 5.5 Funktionstasten und Menüs am Gerätedisplay

Über die Funktionstasten und Menüs des Gerätedisplays können direkt am Autosampler verschiedene Einstellungen vorgenommen und bestimmte Befehle ausgeführt werden.

Allgemeine Informationen zu den Funktionstasten finden Sie im Kapitel 5.5.1 (siehe unten). Informationen zu den einzelnen Menüs finden Sie im Kapitel 5.5.2 ( $\rightarrow$  Seite 76).

#### 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 75).

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



Abb. 29: 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. 30: Funktionstasten

Um	Wählen Sie
das Hauptmenü aufzurufen (→ Seite 78).	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 ( $\rightarrow$ Seite 79).	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 56).	Prime

Wenn der Autosampler in Chromeleon verbunden ist ("connected"), sind keine Eingaben über das Gerätedisplay möglich. Der Drehteller kann jedoch weiterhin über die Taste **Rotate** bewegt werden.

# 5.5.2 Autosampler-Menüs

Abb. 31 zeigt eine Übersicht über die verschiedenen Autosampler-Menüs. Informationen zum allgemeinen Aufbau der Menüs finden Sie auf Seite 77. 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 78).



Abb. 31: Menüstruktur (hier: WPS-3000TPL)

# 5.5.2.1 Aufbau der Menüs

Die einzelnen Menüs sind wie folgt aufgebaut:



Abb. 32: 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

Wählen Sie den gewünschten Menüpunkt über die Pfeiltasten aus—der ausgewählte Menüpunkt ist unterstrichen—und bestätigen Sie die Auswahl mit **Select**. Über **Back** gelangen Sie eine Menüebene zurück.

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 ( $\rightarrow$ <b>Key autorepeat</b> , Seite 81).	^
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 81).	~
zur nächsten Stelle einer Zahl weiterzugehen. Ein eventuell vorhandener Dezimalpunkt 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.	
Hinweis: Abhängig vom ausgewählten Menüpunkt können spezifische Tasten die oben genannten Tasten in der Navigationsleiste ersetzen.	

Wird ein Fehler erkannt, blinken eine oder mehrere Fehlermeldungen auf dem Gerätedisplay. Dann erscheinen folgende Softkeys in der Navigationsleiste:

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 auf.



Abb. 33: Hauptmenü "Main"

Informationen zu den einzelnen Menüs finden Sie in den folgenden Kapiteln:

- Menü Control ( $\rightarrow$  Seite 79)
- Menü Maintenance ( $\rightarrow$  Seite 80)
- Menü Diagnostics ( $\rightarrow$  Seite 81)
- Menü Configuration ( $\rightarrow$  Seite 82)

# 5.5.2.3 Menü Control

Über das Menü **Control** können Sie verschiedene Waschprozeduren ausführen und ein bestimmtes Segment im Drehteller nach vorn holen:

<ul> <li>Control (1.</li> <li>1. Wash</li> <li><u>2. Rotate</u></li> <li>3. Nominal sa</li> <li>4. Sample co</li> </ul>	.4) Imple temp oling		
$\checkmark$	^	Select	Back

Abb. 34: Menü "Control" (hier: WPS-3000TPL)

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
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 ( <b>On</b> oder <b>Off</b> ).	Sample cooling

# 5.5.2.4 Menü Maintenance

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



Abb. 35: Menü "Maintenance"(hier: WPS-3000TPL)

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
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) und können ggf. entsprechende Einstellungen vornehmen oder Befehle ausführen:



Abb. 36: Menü "Diagnostics"

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 ggf. entsprechende Einstellungen vornehmen oder Befehle ausführen:



Abb. 37: Menü "Configuration"

Um	Wählen Sie
<ul> <li>die Einstellungen für das Display und die Funktionstasten festzulegen:</li> <li>Brightness—legt die Helligkeit der Displayanzeige fest (in Prozent).</li> <li>Contrast—legt den Kontrast der Displayanzeige fest (in Prozent).</li> <li>Key sound—legt fest, ob bei Betätigung einer Funktionstaste ein akustisches</li> <li>Signal ertönt (On = ja oder Off = nein).</li> <li>Key autorepeat—legt fest, ob bei längerer Tastenbetätigung der Tastendruck</li> <li>automatisch wiederholt wird, z.B. zur schnellen Änderung eines</li> <li>Einstellwertes (On = ja oder Off = nein).</li> </ul>	Display & soft keys
<ul> <li>Festzulegen, ob und wie eine Leakerkennung erfolgen soll:</li> <li>Enabled—schaltet die Leakerkennung ein. Wird eine Undichtigkeit erkannt, erscheint eine Meldung auf dem Gerätedisplay und es ertönt ein akustischer Alarm.</li> <li>Silent—schaltet die Leakerkennung ein. Wird eine Undichtigkeit erkannt, erscheint nur eine Meldung auf dem Gerätedisplay. Es ertönt kein akustischer Alarm.</li> <li>Disabled—schaltet die Leakerkennung aus.</li> </ul>	Leak sensor mode
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 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
Festlegen der Injektionsmethode	Siehe unten
Synchronisieren von Injektionsbefehl und Kolbenhüben	86
Einschalten der Probenthermostatisierung	86
Ändern der Drehtellerkonfiguration	87
Festlegen der Injektionsparameter und Starten der Injektion	88
Einschalten der Leakerkennung	89
Anpassen von Helligkeit und Kontrast der Displayanzeige 89	
Empfehlungen für den Betrieb	90

# 5.6.1 Festlegen der Injektionsmethode

Der Autosampler unterstützt Full-Loop-, Partial-Loop- und Microliter-Pick-up-Injektionen. Darüber hinaus können Sie in Chromeleon benutzerdefinierte Programme für die Probenvorbereitung und Injektion erstellen. Die Injektionsmethode wird über Chromeleon festgelegt. Nachfolgend ist die Auswahl über den Programm-Assistenten beschrieben. Die genannten Parameter stehen auch im Dialogfenster Commands und auf dem Steuerfenster für den Autosampler zur Verfügung.

- 1. Rufen Sie den Programm-Assistenten auf ( $\rightarrow$  Seite 71).
- 2. Wählen Sie auf der Seite **Sampler Options** unter **Inject Mode** die Injektionsmethode aus: **UserProg**, **Partial**, **FullLoop** oder **ulPickUp**.

**Full-Loop-Injektionen** bieten optimale Reproduzierbarkeit, da die Probenschleife vollständig mit Probe gefüllt wird. Das Injektionsvolumen entspricht dem Probenschleifenvolumen. Der Probenbedarf berechnet sich wie folgt:

FlushVolume + LoopVolume x LoopOverfill

Für alle weiteren Injektionen aus demselben Probengefäß (wenn RinseBetweenReinjections = No): *FlushVolume2* + *LoopVolume x LoopOverfill* 

Informationen zum Injektionsablauf (Verteilung der Flüssigkeiten und Schaltsequenz) finden Sie auf Seite 150.

Bei **Partial-Loop-Injektionen** ist der Probenbedarf geringer, da die Probenschleife nur zum Teil mit Probe gefüllt wird. Außerdem ist das Injektionsvolumen frei wählbar und variabel, ohne dass eine Probenschleife mit einem anderen Volumen installiert werden muss. Der Probenbedarf berechnet sich wie folgt:

#### *FlushVolume* + *InjectionVolume*

Für alle weiteren Injektionen aus demselben Probengefäß: *FlushVolume2* + *InjectionVolume* 

Informationen zum Injektionsablauf (Verteilung der Flüssigkeiten und Schaltsequenz) finden Sie auf Seite 154.

Beim **Mikroliter-Pick-up** wird das Schleifenvolumen vollständig injiziert, enthält aber nur zu einem Teil Probenvolumen. Es wird ein Pfropfen Transportflüssigkeit, gefolgt von dem Probenpfropfen in die Probenschleife transportiert, wiederum gefolgt von Transportflüssigkeit. Daher sollte das TransportVial Elutionsmittel des Startgradienten enthalten. Diese Methode erzeugt ohne Probenverlust höchste Genauigkeit, aber eine etwas geringere Reproduzierbarkeit.

Informationen zum Injektionsablauf (Verteilung der Flüssigkeiten und Schaltsequenz) finden Sie auf Seite 157.

Parameter	Beschreibung
FlushVolume	Legt das Spülvolumen der Nadel fest. Sie können auch ein niedrigeres als das voreingestellte Spülvolumen wählen, mindestens muss dieses jedoch dem Injektionsnadelvolumen entsprechen (Empfehlung: zweifaches Nadelvolumen). Ein niedrigeres Spülvolumen kann allerdings die Messgenauigkeit beeinträchtigen. Für den WPS-3000FC muss das Spülvolumen größer oder gleich der Summe aus dem Nadelvolumen und dem Bridge-Volumen sein (Empfehlung: 2x (Vedelscher und Vedersen))
	(Nadelvolumen + Bridge-Volumen)).
FlushVolume2 (nur bei Full-Loop- und Partial-Loop- Injektionen)	Legt das Spülvolumen der Nadel für weitere Injektionen aus demselben Probengefäß fest. Die Voreinstellung beträgt 2 µL. Voraussetzung ist, dass <b>RinseBetweenReinjections</b> auf <b>No</b> gesetzt ist.
Volume	Legt das Injektionsvolumen fest. Bei Full-Loop-Injektionen sollte <b>Volume</b> dem Volumen der Probenschleife entsprechen.
LoopOverfill (nur bei Full-Loop- Injektionen)	Legt das Volumen fest, das vor der Injektion durch die Probenschleife transportiert wird. Die Voreinstellung ist 2. Dies bedeutet, dass die Probenschleife einmal gespült und dann gefüllt wird.

Informationen zu benutzerdefinierten Injektionsprogrammen finden Sie in Kapitel 5.7.7 ( $\rightarrow$  Seite 96).

3. Die in der nachfolgenden Tabelle genannten Parameter sind für die Verwendung von Wasser als Lösungsmittel optimiert und in Chromeleon voreingestellt. Ändern Sie diese Parameter ggf. bei Verwendung eines anderen Lösungsmittels entsprechend Ihrer Anwendung ab.

Parameter	Beschreibung	Voreinstellung für WPS-3000	
		PL (RS), T(B)PL	TBPL Analytisch / T(B)FC
DispSpeed	Geschwindigkeit, mit der die Probe ausgestoßen wird.	2000 nL/s	20 µL/s
DrawSpeed*	Geschwindigkeit, mit der die Probe aufgezogen wird.	50 nL/s	3 µL/s
DrawDelay	Gibt an, wie lange die Probennadel nach dem Aufziehen der Probe noch im Probenfläschchen bleibt.	5 s	3 s
WashSpeed	Geschwindigkeit der Spritze für den Waschvorgang	4000 nL/s	20 µL/s
WasteSpeed	Geschwindigkeit, mit der die Flüssigkeit aus der Spritze in den Abfall entleert wird.	4000 nL/s	32 µL/s

\* Je nach Gasgehalt der Probe und Injektionsmethode ist auch eine höhere Geschwindigkeit möglich.

# 5.6.2 Synchronisieren von Injektionsbefehl und Kolbenhüben

#### Nur in Verbindung mit einer Pumpe DGP-3600 oder LPG-3400 relevant

Nur wenn Sie den Autosampler im Programm **Server Configuration** mit einer DGP-3600 oder LPG-3400 verbunden haben ( $\rightarrow$  Seite 43), kann der Injektionsbefehl mit den Kolbenhüben der Pumpe synchronisiert werden. Damit wird sichergestellt, dass alle Injektionen zur selben Phase des Pumpenzyklus erfolgen. Auf diese Weise wird bei Gradientenanwendungen die Retentionszeitpräzision deutlich verbessert.

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

- Setzen Sie SyncWithPump auf Off, wenn Sie ohne Synchronisierung arbeiten möchten.
- 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 gewünschten 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 können Sie über Chromeleon oder am Gerätedisplay ein- und ausschalten. Ebenso können Sie dort die gewünschte Temperatur festlegen.

#### Über Chromeleon

- Öffnen Sie in Chromeleon das Dialogfenster Commands f
  ür den Autosampler (→ Seite 68).
- 2. Wählen Sie **Temperature** und **Nominal** und geben Sie die gewünschte Temperatur ein. Durch die Eingabe einer Temperatur wird **TempCtrl** auf **On** gesetzt, falls noch nicht geschehen.



Abb. 38: Einschalten der Probenthermostatisierung

Setzen Sie **TempCtrl** auf **Off**, wenn Sie für eine bestimmte Anwendung ohne Probenthermostatisierung arbeiten möchten.

Wenn Sie die Thermostatisierung 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.

### Am Gerätedisplay

- 1. Rufen Sie am Gerätedisplay das Menü Control auf ( $\rightarrow$  Seite 79).
- 2. Wählen Sie **Nominal sample temperature** und geben Sie die gewünschte Temperatur ein.
- 3. Wählen Sie Sample Cooling und On, um die Probenthermostatisierung einzuschalten.

Setzen Sie Sample Cooling auf Off, wenn Sie ohne Thermostatisierung 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 gewünschte Änderung im Programm Server Configuration auf der Seite Segments / Pump Link vor (→ Seite 43).
- Wenn Sie die Standardeinstellung nicht ändern möchten, können Sie 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. Um dies zu vermeiden, können Sie den automatischen Tray-Test aktivieren ( $\rightarrow$  Seite 95).

# 5.6.5 Festlegen der Injektionsparameter und Starten der Injektion

In Chromeleon können Sie die Injektion manuell oder mit Hilfe eines Programms durchführen. In jedem Fall sendet der Autosampler ein Signal an Chromeleon, wenn die Injektion erfolgt ist.

#### Manuelle Injektion

■ Hinweis: Vergewissern Sie sich zunächst, dass die gewünschte Injektionsmethode ausgewählt ist (→ Seite 83).

Wählen Sie eine der folgenden Alternativen, um die Injektionsparameter einzugeben und die Injektion zu starten:

• Dialogfenster Commands

Wählen Sie **Inject**, legen Sie unter **Position** die Probenposition und unter **Volume** das Injektionsvolumen fest. Klicken Sie dann auf die Schaltfläche **Execute**, um die Injektion zu starten.

• Steuerfenster für den Autosampler

Legen Sie im Abschnitt **Inject** die Probenposition und das Injektionsvolumen fest und starten Sie die Injektion über die Schaltfläche **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 und klicken Sie **OK**, um die Injektion zu starten.

## Programmgesteuerte Injektion

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 73).

# 5.6.6 Einschalten der Leakerkennung

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

#### Einschalten der Leakerkennung über Chromeleon

- Öffnen Sie in Chromeleon das Dialogfenster Commands f
  ür den Autosampler (→ Seite 68).
- 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** ( $\rightarrow$  Seite 74).
- Rufen Sie das Menü Configuration auf und wählen Sie den Punkt Leak sensor mode (→ Seite 82).
- 3. Wählen Sie **Enabled** oder **Silent** (siehe oben). Über **Disabled** wird die Leakerkennung ausgeschaltet.

## 5.6.7 Anpassen von Helligkeit und Kontrast der Displayanzeige

Die Helligkeit der Hintergrundbeleuchtung und den Kontrast der Displayanzeige können Sie, falls erforderlich, 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 und/oder unter Contrast den Wert für den Kontrast der Displayanzeige
- Rufen Sie am Gerätedisplay das Menü Configuration und dann Display & soft keys auf (→ Seite 81). Ändern Sie unter Brightness den Wert für die Helligkeit der Hintergrundbeleuchtung und/oder unter Contrast den Wert für den Kontrast der Displayanzeige.

# 5.6.8 Empfehlungen für den Betrieb

Es wird empfohlen, folgende Aktionen in regelmäßigen Abständen, besonders nach längeren Stillstandszeiten, durchzuführen:

- Prüfen Sie, ob sich Luftblasen in der Spritze befinden und spülen Sie die Spritze, falls erforderlich (→ Seite 56).
- Führen Sie einen Waschzyklus für die Pufferschleife durch.
   Führen Sie dazu den Befehl Wash in Chromeleon aus oder wählen Sie am Gerätedisplay die Funktionstaste Wash (→ Seite 74), oder wählen Sie Wash im Menü Control (→ Seite 79). Zum Waschen der Pufferschleife sollte mindestens das doppelte Volumen der Pufferschleife verwendet werden. Dazu muss die Spritze gegebenenfalls öfters aufgezogen werden.
- Prüfen Sie die Probennadel in regelmäßigen Abständen auf Abnutzung und Beschädigung. Dies ist besonders wichtig bei Probennadeln aus Fused Silica oder PEEK.

# 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
Definieren der Nadelhöhe und Vorstecher-Einstechtiefe (Sample Height, Trans Liquid Height)	93
Bewegen des Drehtellers vor dem Aufziehen der Probe (Tray Shake)	94
Operational Qualification und Performance Qualification	95
Automatischer Tray Test	95
Automatische Probenfläschchenerkennung	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 Hilfe zu Chromeleon.

# 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 gewünschten Grenzwerte fest. Eine vollständige Liste der für den Autosampler verfügbaren Befehle und Zähler finden Sie in der Hilfe zu Chromeleon. Um die Informationen für die Predictive Performance aktuell zu halten, werden folgende Maßnahmen empfohlen ( $\rightarrow$  Tabelle).

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

Damit werden die entsprechenden Zähler zurückgesetzt. Außerdem wird das Datum eingetragen, an dem die Maßnahme erfolgt ist.

#### Steuerfenster

Auf dem Steuerfenster für den Autosampler ( $\rightarrow$  Seite 69) 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 gegebenenfalls 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, und es ist nicht möglich, einen Batch zu starten.

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

# 5.7.2 Definieren der Nadelhöhe und Vorstecher-Einstechtiefe

Der Parameter **SampleHeight** beschreibt den Abstand (in mm) zwischen dem äußeren Boden des Probengefäßes (Mikrotiterplatte, Deep-Well-Plate oder Probenfläschchen) und der Nadelspitze und definiert so, wie tief die Probennadel beim Ansaugen der Probe in das Probengefäß eintaucht. Die Nadelhöhe für die Reagens- und Transportfläschchen wird analog über den Parameter **TransLiquidHeight** eingestellt. Beachten Sie für die Einstellung der Nadelhöhe auch die Informationen auf Seite 94.

Der Parameter **PunctureDepth** definiert, wie tief (in mm) der Vorstecher in das Probengefäß (Mikrotiterplatte, Deep-Well-Plate oder Probenfläschchen) einsticht. Dies wird vom oberen Rand des Probengefäßes gemessen. Die Einstechtiefe für die Reagensund Transportfläschchen wird analog über den Parameter **TransVialPunctureDepth** eingestellt. Beachten Sie für die Einstellung der Einstechtiefe auch die Informationen auf Seite 94.





Nr.	Beschreibung
1	Nadelpaar (Vorstecher und Probennadel)
2	Einstechtiefe des Vorstechers (PunctureDepth, TransVialPunctureDepth)
3	Probengefäß
4	Höhe der Probennadel über dem inneren Boden des Probengefäßes (SampleHeight, TransLiquidHeight)
5	Ausgleichswert für die Bodenhöhe (SampleHeightOffset_XX)
6	Probenhalter
7	Drehteller

Beachten Sie bei der Einstellung der Nadelhöhe folgende Hinweise:

- Setzen Sie den Parameter **SampleHeight** *nicht* auf '**0**', wenn Sie ein Probengefäß installiert haben. Bei dieser Einstellung berührt die Probennadel den Boden des Probengefäßes und kann beschädigt werden.
- Wenn Sie aus einem Probengefäß nur *ein* Mal injizieren möchten, sollten Sie **SampleHeight** so wählen, dass die Probennadel nur so tief wie nötig in das Probengefäß eintaucht. Damit wird eine Verschleppung zwischen den Proben weitgehend vermieden.
- Wenn Sie *mehrfach* Probe aus einem Probengefäß aufziehen möchten, sollten Sie **SampleHeight** so wählen, dass die Probennadel so tief wie möglich, also bis knapp über den Boden, in das Probengefäß eintaucht. Damit wird sichergestellt, dass die Probe möglichst vollständig aufgezogen werden kann.
- Um zu verhindern, dass die Probennadel den Boden des Probengefäßes berührt, wenn **SampleHeight** auf 0 mm gesetzt ist, steht der Parameter **SampleHeightOffset\_XX** zur Verfügung. Geben Sie hier die Bodenhöhe des installierten Probengefäßtyps ein, damit diese mit berücksichtigt wird. Standardmäßig ist der Wert für **SampleHeightOffset\_XX** auf null gesetzt.
- Bei sehr kleinen Einstellungen für **TransLiquidHeight** müssen Sie unter Umständen die **TransVialPunctureDepth** erhöhen, um den Boden des Transportfläschchens erreichen zu können.

Beachten Sie bei der Einstellung der Einstechtiefe für den Vorstecher folgende Hinweise:

- Wählen Sie die Einstechtiefe so, dass das Septum des Probengefäßes durchstochen wird, der Vorstecher aber noch nicht in die Probe eintaucht.
- Wird die Einstechtiefe zu niedrig gewählt, wird das Septum des Probengefäßes nicht durchstochen. Die Probennadel gelangt nicht in das Probengefäß und kann beschädigt werden.
- Für die meisten HPLC-Probenfläschchen, ist eine Einstechtiefe von 8 mm ausreichend, damit das Septum des Probenfläschchens durchstochen wird.

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

Zur Re-Homogenisierung der Probe, speziell nach längeren Standzeiten, unterstützt Chromeleon den "Tray Shake". Durch schnelle Drehtellerbewegungen wird die Probe dabei vor dem Aufziehen gleichmäßig durchmischt. Peakflächen-Änderungen aufgrund einer lokalen Änderung der Konzentration im Probengefäß werden so verhindert.

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.

# 5.7.4 Operational Qualification und Performance Qualification

Sie können die Leistungsfähigkeit des Systems mit Hilfe der Operational und Performance Qualification kontrollieren und dokumentieren. Alle erforderlichen Materialien sowie eine detaillierte Anleitung zur Durchführung sind auf Anfrage erhältlich.

# 5.7.5 Automatischer Tray-Test

Diese Funktion ist standardmäßig deaktiviert. Sie kann jedoch unter Chromeleon eingeschaltet werden. Setzen Sie dazu das Property EnableTravTest auf Enabled.

Der automatische Tray-Test überprüft, 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 Trav-Test wird vor jeder ersten Injektion aus einem Probenhalter nach einem manuellen Eingriff einmal pro Probenhalter ausgeführt.

## 5.7.6 Automatische Probengefäßerkennung

Ein Sensor im Niederhalter dient zur automatischen Probenfläschchen- bzw. Mikrotiterplatten- und Deep-Well-Plate-Erkennung.



Abb. 40: Niederhalter

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

**Hinweis:** Für 24-Deep-Well-Plates ist ein Adapter am Niederhalter erforderlich (Best.-Nr. 6820.2402). Ist der Adapter installiert, darf der automatische Tray-Test nicht aktiviert sein ( $\rightarrow$  Kapitel 5.7.5).

# 5.7.7 Benutzerdefinierte Programme

### 5.7.7.1 Benutzerdefinierte Programme erstellen

Mit dem Programm-Assistenten von Chromeleon können Sie neben Programmen zur Probenanalyse Chromeleon auch benutzerdefinierte Programme erstellen, mit denen Sie beispielsweise die einzelnen Schritte zur Probenvorbereitung genau festlegen können.

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

ogram Wizard: Sampler Option	S		
Inject Mode:	UserProg	•	
	UserProg	1	
Transport Vials (µIPickUp):	FullLoop		To: RA1
Transport Vial Capacity:	Unlimited		[099999]
Transport Liquid Height:	5,000		[0,00030,000 mm]
Transport Vial Puncture Depth:	5,000		[0,00011,000 mm]
	2 400		70 400 40000 000 II
riusri volume (ruiiLoop/raniai).	2,400		[2,40010000,000 µ]
Flush Volume 2:	2,000		[0,00010000,000 µl]
Loop Overfill:	2,000		[1,00010,000]

Abb. 41: Benutzerdefiniertes Programm

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

Das benutzerdefinierte Programm wird dann mit dem Befehl **Inject** des Programms zur Probenanalyse gestartet. Der chromatographische Lauf wird angehalten, bis über das benutzerdefinierte Programm der Befehl UDPInjectMarker gegeben wird.

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

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

# 5.7.7.2 Eingabe relativer Probenpositionen

Chromeleon unterstützt in benutzerdefinierten Programmen auch die Angabe von relativen Probenpositionen, beispielsweise für das Probenfläschchen, dass zur Probenvorbereitung verwendet werden soll (PrepVial).

Dabei kann sich die so angegebene Position in einer anderen Reihe auf dem Probenhalter, in einem anderen Drehtellersegment, oder eine bestimmte Anzahl von Positionen von einem anderen Probenfläschchen entfernt befinden.

Informationen darüber, wie Sie relative Probenpositionen in einem Programm eingeben müssen und was Sie darüber hinaus beachten sollten, finden Sie in der Hilfe zu Chromeleon.

# 5.8 Außerbetriebnahme des Autosamplers

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

- Spülen Sie den Autosampler von Lösungsmitteln frei. Wird ein Puffer als Teil der mobilen Phase eingesetzt, spülen Sie das System vor der Außerbetriebnahme einige Male mit einem Gemisch aus Methanol und Wasser (50:50), um eine Aufkonzentrierung von Salzen im System zu vermeiden.
- Füllen Sie den Autosampler mit Methanol oder einem ähnlichen Alkohol wie Isopropanol oder Ethanol. Sind die im Autosampler enthaltenen Lösungsmittel nicht wasserlöslich, müssen diese schrittweise ersetzt werden.
- Wenn Sie den Autosampler transportieren oder versenden möchten, müssen Sie
  - die beweglichen Teile des Drehtellers (Probenhalter, Probenfläschchen, Mikrotiterplatten) und gegebenenfalls die Flasche für die Waschflüssigkeit entnehmen.
  - den Nadelarm und den Drehtellerdeckel wie in der Verpackungsvorschrift beschrieben über das entsprechende Schaumteil sichern.
  - ▲ Vorsicht: Dies ist ein Muss! Um eine Beschädigung der Probennadel und des Waschports zu vermeiden, fahren Sie die Probennadel und den Nadelarm zuerst in die korrekte Position (Nadel rechts und Nadel nicht im Waschport). Drücken Sie dazu 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 und beachten Sie die Verpackungsvorschrift. Erfolgt der Versand des Gerätes 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 und auch auf Anfrage erhältlich.

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

#### Standby-Programm

Ein Standby-Programm versetzt die HPLC-Anlage in den Bereitschaftszustand. Dadurch kann die jeweilige Applikation schnell wieder aktiviert werden. Ein Standby-Programm umfasst im Wesentlichen die folgenden Punkte:

- Der Fluss wird am Programmende automatisch verringert.
- Die Temperatur aller temperaturgesteuerten Systemmodule wird heruntergefahren.

#### Shutdown-Programm

Ein Shutdown-Programm fährt die HPLC-Anlage automatisch herunter. Es umfasst im Wesentlichen die folgenden Punkte:

- Der Fluss wird am Programmende automatisch gestoppt.
- Bestimmte Systemfunktionen werden abgeschaltet (z.B. Detektorlampen, Temperaturregelung der temperaturgesteuerten Systemmodule).

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

Wählen Sie eine der folgenden Alternativen:

- Wählen Sie die erforderlichen Befehle und Parameter im Dialogfenster Commands aus (→ Seite 68).
- Automatisieren Sie die Außerbetriebnahme, indem Sie ein entsprechendes Programm erstellen und ablaufen lassen ( $\rightarrow$  Seite 71).
  - Hinweis: Für den WPS-3000TBPL Analytisch in einem analytischen UltiMate 3000-System steht Ihnen in Chromeleon zusätzlich der SmartShutdown-Assistent zur Verfügung. Mit Hilfe dieses 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. Weitere Informationen finden Sie in der Hilfe zu Chromeleon

# 5.9 Wartung und Wartungsintervalle

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

Die unten genannten Wartungsarbeiten sollten Sie in regelmäßigen Intervallen durchführen, um die optimale Funktionsfähigkeit und maximale Verfügbarkeit Ihres Autosamplers sicherzustellen. Dabei soll Ihnen die untenstehende Tabelle als Orientierungshilfe dienen, 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	Überprüfen Sie vor Betriebsbeginn die fluidischen Verbindungen auf mögliche Luftblasen und entgasen Sie die Waschflüssigkeit.
	Überprüfen Sie die Fluidik auf mögliche Undichtigkeiten.
	Überprüfen Sie die fluidischen Verbindungen auf mögliche Salzablagerungen.
	Wenn Sie mit Pufferlösungen arbeiten, sollten Sie das System nach Abschluss der Arbeiten gründlich mit einer Flüssigkeit spülen, die keine Puffer/Salze enthält.
Regelmäßig	Überprüfen Sie die Schlauchverbindungen auf mögliche Beschädigungen (Knicke, Risse, Schnitte, Blockierung).
	Überprüfen Sie den Drehteller auf Rückstande und entfernen Sie diese ggf. mit einem Tuch oder Papier. Entnehmen Sie dazu die Probenhalter und, falls erforderlich, die 5-Positionen Halter. Diese Überprüfung ist speziell für den Autosampler mit Probenthermostatisierung wichtig, da ein sauberer Drehteller die Kühlleistung verbessert.
	Überprüfen Sie die Schläuche, die zur Ableitung von Flüssigkeit aus dem Innenraum an die beiden Abläufe rechts unterhalb des Autosamplers angeschlossen sind (→ Seite 56). Die Schläuche dürfen nicht abgeknickt sein und an keiner Stelle höher als der Anschluss-Stutzen liegen. Entleeren Sie den Abfallbehälter, wenn erforderlich.
Häufigkeit	Was ist zu tun
------------	--
Jährlich	<ul> <li>Überprüfen Sie folgende Komponenten und tauschen Sie diese ggf. aus:</li> <li>Verbindungen an den Schaltventilen (→ Seiten 55 bzw. 137)</li> <li>Verbindungen an der Spritze (→ Seite 55)</li> <li>Probennadel (→ Seite 121)</li> <li>Vorstecher (→ Seite 123)</li> <li>Rotordichtung und Stator (→ Seite 125)</li> </ul>
	Lassen Sie den Thermo Fisher Scientific-Kundendienst eine regelmäßige Überprüfung auf Verschmutzung, Abnützung, usw. in Abständen von ca. 1 Jahr durchführen

■ Hinweis: Unter Chromeleon stehen Funktionen zur Lebensdauerprognose von Verschleißteilen zur Verfügung (→ Seite 91).

# 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:

#### Statusanzeigen

Die Statusanzeigen 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 21).

#### Fehlermeldungen

Wird während des Betriebs des Autosamplers ein Fehler erkannt, erscheint eine entsprechende Fehlermeldung auf dem Gerätedisplay. Im Kapitel 6.2 ( $\rightarrow$  Meldungen, Seite 104) finden Sie zu jeder Meldung eine kurze Beschreibung der möglichen Ursachen sowie Vorschläge für Abhilfemaßnahmen. Wird der Autosampler unter Chromeleon betrieben, wird der Fehler auch im Chromeleon Audit Trail protokolliert.

**Hinweis:** Hinweise zu Störungen, die beim Betrieb des UltiMate 3000 RSLCnano-Systems auftreten können, finden Sie im Handbuch für das NCS Nano-Cap-System.

> Hinweise zu Störungen, die beim Betrieb des UltiMate 3000-Systems mit Flow Manager für Anwendungen in der Nano- oder Kapillar-HPLC auftreten können, finden Sie im UltiMate 3000 Proteomics MDLC-Systemhandbuch.

> Hinweise zu Störungen, die beim Betrieb des UltiMate 3000-Systems für Anwendungen in der analytischen HPLC auftreten können, finden Sie in den Handbüchern zu den anderen Modulen des Systems.

Wenn Sie den Fehler nicht mit Ihren Mitteln beheben können, wenden Sie sich bitte an den Thermo Fisher Scientific-Kundendienst.

# 6.2 Meldungen auf dem Gerätedisplay

Tritt während des Betriebs des Autosamplers ein Fehler auf, leuchtet die LED **Status** auf der Gerätevorderseite rot und es erscheinen eine oder mehrere Meldungen auf dem Autosampler-Display. In der Navigationsleiste erscheinen dann 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. Beachten Sie, dass einige der unten genannten Meldungen für die verschiedenen Antriebe identisch sind. 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 Meldungen auf dem Gerätedisplay auch über den Chromeleon-Befehl **ClearDisplayError** gelöscht werden.

Die nachfolgende Tabelle listet Meldungen auf, die beim Betrieb des Autosamplers auftreten können, und nennt entsprechende Abhilfemaßnahmen. Zusätzlich zu den genannten Meldungen können noch weitere Meldungen erscheinen. Wenden Sie sich in diesem Fall bitte mit dem genauen Wortlaut der Meldung an den Thermo Fisher Scientific-Kundendienst.

Meldung	Abhilfe
24V power failure.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): End point not reached.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite 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 Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Home position not found.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Home sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Needle sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.

Meldung	Abhilfe
(Antrieb): No communication.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Not initialized.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Not ready.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
(Antrieb): Obstruction detected.	<ul> <li>Es befindet sich ein Hindernis am Niederhalter oder die Probennadel wird beim Eintauchen in das Probengefäß blockiert. Entfernen Sie das Hindernis.</li> <li>Die Probennadel sticht auf den Rand des Probengefäßes, nicht in das Septum. Überprüfen Sie, ob die Angaben auf der Registerkarte Segments / Pump Link in Chromeleon (→ Seite 43) mit der tatsächlichen Bestückung des Drehtellers übereinstimmt.</li> </ul>
(Antrieb): Stop condition at start.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
A self test is required.	Es muss ein Selbsttest durchgeführt werden. Der Autosampler funktioniert erst wieder, wenn der Selbsttest erfolgreich durchgeführt wurde. 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 die entsprechende Meldung.
Buffer loop is too full.	Die Pufferschleife enthält zu viel Flüssigkeit. Leeren Sie die Pufferschleife über die Funktionstaste <b>Wash</b> ( $\rightarrow$ Seite 74) oder den Befehl <b>Wash</b> im Menü <b>Control</b> ( $\rightarrow$ Seite 79).
Carousel temperature sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Command input overflow.	Führen Sie ein Firmware- oder Treiberupdate durch.
Command queue overflow.	Das benutzerdefinierte Programm ist zu lang. Kürzen Sie das Programm.
Command syntax error.	Führen Sie ein Firmware- oder Treiberupdate durch.
Command was aborted.	Der Befehl wurde durch den Anwender 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 aus und wieder ein oder führen Sie einen Selbsttest durch.
Device is off.	Der Autosampler ist ausgeschaltet (Standby). Schalten Sie den Autosampler über die Standby-Taste auf der Gerätevorderseite oder in Chromeleon ein.
DVC position error.	Vergewissern Sie sich, dass das Injektionsventil nicht blockiert ist.

Meldung	Abhilfe
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 Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Illegal command parameter.	Führen Sie ein Firmware- oder Treiberupdate durch.
Illegal syringe speed.	Die angegebene Spritzengeschwindigkeit ist nicht zulässig. Ändern Sie die Spritzengeschwindigkeit.
Illegal value.	Führen Sie ein Firmware- oder Treiberupdate durch.
Inject volume must be equal to loop volume.	Das Injektionsvolumen entspricht nicht dem Volumen der Probenschleife. (Die Meldung tritt nur im Full-Loop-Injektionsmodus auf.) Ändern Sie das Injektionsvolumen.
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 für die Mischoperation angegebene Ziel ist nicht gültig. Geben Sie ein gültiges Ziel ein (Chromeleon-Befehle UdpDraw > From bzw. UdpDispense > To).
Leak detected.	Der Leaksensor des Autosamplers hat eine Undichtigkeit erkannt. Prüfen Sie, an welcher Stelle Flüssigkeit austritt und beheben Sie die Ursache für die Undichtigkeit. Trocknen Sie den Leaksensor und die Auffangwanne $(\rightarrow \text{Kap. 7.2, Seite 112}).$
Motor switching valve malfunction.	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 Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Needle home position not found.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Needle home sensor malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Needle malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Out of transport liquid.	Es ist keine Flüssigkeit mehr in den Transportfläschchen vorhanden. Befüllen Sie die Transportfläschchen neu und setzen Sie in Chromeleon den Zähler <b>TransportVialUsecount</b> auf null zurück.

Meldung	Abhilfe
Property is protected.	Sie haben versucht, eine geschützte Variable zu ändern. Einige Variablen sind geschützt und können vom Anwender nicht geändert werden.
Property is read only.	Sie haben versucht, eine Lesevariable zu ändern. Einige Variablen sind Lesevariablen und können vom Anwender nicht geändert werden.
Reagent vial not found.	An der angegebenen Position befindet sich kein Reagensfläschchen. Stellen Sie ein Reagensfläschchen an die angegebene 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.	<ul> <li>Die Probennadel taucht nicht tief genug in das Probengefäß ein und kann keine Flüssigkeit aufziehen. Erhöhen Sie, wenn möglich, die Einstechtiefe in Chromeleon</li> <li>(= PunctureDepth für Probengefäße bzw. TransVialPunctureDepth für Transport- und Reagensfläschchen).</li> <li>Befüllen Sie alternativ das Probengefäß neu oder ändern Sie die Einstellung für die Nadelhöhe (→ Seite 93). Evtl. ist der Wert SampleHeightOffset_XX nicht auf null gesetzt.</li> </ul>
Syntax error.	Führen Sie ein Firmware- oder Treiberupdate durch.
Syringe is not empty.	Die Spritze ist zu Beginn einer Injektion nicht leer. 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 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.	Überprüfen Sie das Spritzenventil.
Temperature difference too large. Probable coolant circuit malfunction.	Schließen Sie ggf. den Drehteller-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 des Kommandos beendet ist. Führen Sie das gewünschte Kommando erneut aus.
The volume is too large.	Es soll mehr Volumen aufgezogen werden, als die Spritze oder Pufferschleife fassen kann. Ändern Sie das Volumen.
Transport liquid vial not found.	An der angegebenen Position befindet sich kein Transportfläschchen. Stellen Sie ein Transportfläschchen an die angegebene Position oder geben Sie eine andere Position an.
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 Sie die Positionsangabe. Prüfen und korrigieren Sie ggf. die Segmenteinstellungen (→ Registerkarte Segments / Pump Link, Seite ).
USB interface is not ready.	Überprüfen Sie die USB-Verbindung. Ersetzen Sie ggf. ein fehlerhaftes USB-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 der Test zur Probengefäßerkennung ( $\rightarrow$ Seite 95) ausgeschaltet ist und Ihre Probengefäße nicht die erforderliche Mindesthöhe haben ( $\rightarrow$ Seite 24). Thermo Fisher Scientific empfiehlt die Verwendung der auf Seite 24 genannten Probengefäße.
	Es ist ein falscher Segmenttyp definiert. Prüfen und korrigieren Sie ggf. die Segmenteinstellungen (→ Registerkarte Segments / Pump Link, Seite 43).
Vial pusher malfunction.	Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite aus und wieder ein oder führen Sie einen Selbsttest durch.
Wrong tray type detected.	<ul> <li>Bei der automatischen Probenhaltererkennung wurde festgestellt, dass die Angaben auf der Registerkarte Segments / Pump Link nicht mit der tatsächlichen Bestückung des Drehtellers übereinstimmen. Prüfen und korrigieren Sie ggf. die Angaben auf der Registerkarte Segments / Pump Link in Chromeleon (→ Seite 43).</li> <li>Prüfen Sie, ob Ihre Probengefäße die erforderliche Mindesthöhe (→ Seite 24) haben. Thermo Fisher Scientific empfiehlt die Verwendung der auf Seite 24 genannten Probengefäße.</li> </ul>

Wenn der Autosampler unter Chromeleon betrieben wird und keine Kommunikation zwischen dem Autosampler und Chromeleon aufgebaut werden kann, können entsprechende Fehlermeldungen im Chromeleon Audit Trail erscheinen. (Die genannte Nummer 1610103 ist in diesem Fall die Seriennummer des Geräts.)

Meldung	Abhilfe	
WPS-3000PL@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-3000PL@ 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-3000PL@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 gegebenenfalls den in der Meldung angegebenen Autosampler aus der Konfiguration oder wählen in der Serverkonfiguration einen anderen Autosampler aus.	
Error reading from WPS- 3000PL @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 (→ Seite 34). Ersetzen Sie ein evtl. fehlerhaftes USB-Kabel oder Hub.	
Error reading from WPS- 3000PL @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.	

# 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.
- Der Autosampler darf nicht mit offenem Gehäuse betrieben werden.
- Verwenden Sie ausschließlich die von Thermo Fisher Scientific für das Gerät autorisierten Original-Ersatzteile.
- Falls der Autosampler 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 und auf Anfrage erhältlich.

Hinweise zur Außerbetriebnahme des Autosamplers erhalten Sie auf Seite 97.

## 7.2 Leaksensor

Der Leaksensor spricht an, wenn er Flüssigkeit ausgesetzt ist. Beseitigen Sie die Ursache für die Undichtigkeit und trocknen Sie den Leaksensor (tragen Sie dabei geeignete Schutzkleidung):

- 1. Schalten Sie den Autosampler aus.
- 2. Überprüfen Sie den Autosampler auf eventuelle Undichtigkeit. Ziehen Sie undichte Schlauch- und Kapillarverbindungen nach oder tauschen Sie diese aus.
- 3. Entfernen Sie die Flüssigkeit und trocknen Sie den Leaksensor. Saugen Sie dazu die Flüssigkeit, die am unteren Ende des Leaksensors in der Auffangwanne steht, mit einem Tuch auf.



Achten Sie darauf, die Sensoren nicht zu verbiegen oder zu beschädigen.



Abb. 42: Trocknen des Leaksensors

- 4. Der Sensor braucht einige Minuten Zeit, sich auf die Umgebungstemperatur einzustellen.
- 5. Schalten Sie den Autosampler an.
- 6. Wird nach dem Einschalten des Autosamplers kein Fehler gemeldet, kann der Betrieb wieder aufgenommen werden.
- Hinweis: Die LED Status auf der Gerätevorderseite bleibt rot, bis der Sensor trocken ist. Wenn eine Meldung auf dem Gerätedisplay erschienen ist, können Sie diese über die Taste Clear auf der Navigationsleiste löschen (→ Seite 78).

Beschreibung	BestNr.
WPS-3000(T)PL (RS) Nano/Cap, WPS-3000TBPL Nano/Cap	6822 0001
WPS-3000TBPL Analytisch	(822.0001
$100 \ \mu\text{L-Spritze}$ WPS-3000T(B)FC	6822.0002
250 µL-Spritze	6822.0003
Informationen zu optional erhältlichen Spritzen finden Sie im Kapitel 10.2 (→ Seite 171).	

## 7.3 Wechseln der Spritze

- 1. Geben Sie das Kommando **InitiateChangeSyringe** über Chromeleon— *oder* rufen Sie am Gerät das Menü **Maintenance** auf und wählen Sie dann **Syringe** und **Down**.
- 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)
  und entfernen Sie die Schraube dann per Hand. Heben Sie die Schraube f
  ür den
  Einbau der neuen Spritze auf.



Abb. 43: Lösen der Spritze am Spritzenstößel

4. Schrauben Sie die Spritze am Spritzenventil ab.



Abb. 44: Lösen der Spritze am Spritzenventil

5. Vergewissern Sie sich, dass beim Ausbau der alten Spritze die Dichtung im Ventil verblieben ist. (Dies können Sie leicht feststellen, wenn Sie mit einem Spiegel in den unteren Anschlussport des Ventils schauen.)

Falls die Dichtung zusammen mit der alten Spritze entfernt oder beschädigt wurde, installieren Sie bitte eine neue Dichtung, ehe Sie die neue Spritze anschließen. Die Dichtung ist im Lieferumfang der jeweiligen Ersatzspritze enthalten; sie ist jedoch 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. Befüllen Sie die neue Spritze mit Hilfe einer Plastikspritze oder gehen Sie wie unten beschrieben vor:
  - a) Befüllen 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 und schrauben Sie die Spritze dann am Spritzenventil fest.



Abb. 45: Flüssigkeitsfilm am Spritzeneingang

- 10. Richten Sie den Spritzenstößel aus und schrauben Sie die Befestigungsschraube mit der Hand ein. Ziehen Sie die Schraube dann mit dem Maulschlüssel etwas mehr als handfest an.
- 11. Klappen Sie den Frontdeckel nach unten.

- 12. Geben Sie das Kommando **TerminateChangeSyringe** über Chromeleon— *oder* rufen Sie am Gerät das Menü **Maintenance** auf und wählen Sie dann **Syringe** und **Up**.
  - 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 wie im Kapitel 4.6 (→ Seite 56) beschrieben. Denken Sie daran, dass Luft in der Spritze die Messergebnisse stark beeinflusst. Vergewissern Sie sich daher, dass die Spritze luftfrei ist, ehe Sie mit Ihrer Analyse beginnen.
  - **Hinweise:** Wenn die neue Spritze ein anderes Volumen hat als die Spritze, die Sie ausgebaut haben, sollten Sie sich vergewissern, dass
    - in der Konfiguration des Autosamplers auf der Seite Options die Spritzengröße korrekt angegeben ist (→ Seite 44). Ändern Sie die Einstellung gegebenenfalls und übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Dazu steht Ihnen auf der Seite General die Schaltfläche Send configuration to module zur Verfügung (→ Seite 41).
    - auch die passenden Puffer- und Probenschleifen installiert sind (→ Seite 16).

Aktualisieren Sie nach einem Spritzenwechsel die entsprechenden Serviceinformationen in Chromeleon über den Befehl SyringeChanged.

## 7.4 Wechseln der Pufferschleife

Beschreibung	BestNr.
WPS-3000(T)PL (RS) Nano/Cap 50 μL-Pufferschleife	6820.0019
<i>WPS-3000TBPL Nano/Cap</i> 50 μL-Pufferschleife	6821.0019
WPS-3000TFC 500 μL-Pufferschleife	6820.0020
WPS-3000TBPL Analytisch und WPS-3000TBFC 500 µL-Pufferschleife	6821.0020
Informationen zu optional erhältlichen Pufferschleifen finden Sie im Kapitel 10.2 ( $\rightarrow$ Seite 171).	

- 1. Klappen Sie den Frontdeckel nach oben auf.
- 2. Schrauben Sie die Pufferschleife am Spritzenventil ab.



Abb. 46: Pufferschleife am Spritzenventil

- 3. Schrauben Sie die Pufferschleife am Injektionsventil ab (Port 1).
- 4. Schließen Sie die neue Pufferschleife folgendermaßen an, um zu verhindern, dass Luft in das System gelangt:
  - a) Führen Sie über Wash im Menü Control (→ Seite 79) oder in Chromeleon (Wash und WashVolume) einen Waschvorgang durch. Zum Waschen der Pufferschleife sollte mindestens das doppelte Volumen der Pufferschleife verwendet werden. Schließen Sie die neue Pufferschleife an, während der Waschvorgang ausgeführt wird.
  - b) Schrauben Sie die neue Pufferschleife am Spritzenventil fest, sobald die Waschflüssigkeit am Spritzenventil austritt. Dies verhindert, dass beim Einbau der Pufferschleife Luft in die Schleife eindringt.
  - c) Schrauben Sie die Pufferschleife am Injektionsventil fest, sobald die Waschflüssigkeit am freien Ende der Pufferschleife austritt.
- 5. Beachten Sie die Hinweise auf der nächsten Seite.

**I** Hinweise: Wenn die neue Pufferschleife ein anderes Volumen hat als die Pufferschleife, die Sie ausgebaut haben, sollten Sie sich vergewissern, dass

- in der Konfiguration des Autosamplers auf der Seite Options das Volumen der Pufferschleife korrekt angegeben ist (→ Seite 44). Ändern Sie die Einstellung gegebenenfalls und übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Dazu steht Ihnen auf der Seite General die Schaltfläche Send configuration to module zur Verfügung (→ Seite 41).
- auch eine geeignete Probenschleife und Spritze installiert sind (→ Seite 16).

# 7.5 Nadelpaar

Das Nadelpaar besteht aus Vorstecher und Probennadel.



Nadelpaar bestehend aus Vorstecher und Probennadel.

Abb. 47: Nadelpaar

#### • Vorstecher

Der Vorstecher ist eine hohle Stahlnadel, in der sich die Probennadel befindet. Er durchsticht das Septum/die Versiegelung des Probengefäßes, ehe die Probennadel in das Gefäß eintaucht. Informationen zum Wechseln des Vorstechers finden Sie im Kapitel 7.5.4 ( $\rightarrow$  Seite 123).

#### • Probennadel

Die Probennadel befindet sich im Vorstecher. Sie nimmt die Probe auf und transportiert sie in die Probenschleife. Informationen zum Wechseln der Probennadel sowie zu optional verwendbaren Kapillaren finden Sie im Kapitel 7.5.3 ( $\rightarrow$  Seite 121).

## 7.5.1 Überprüfen der Nadelpaar-Position

Gehen Sie folgendermaßen vor, wenn Sie prüfen möchten, ob das Nadelpaar korrekt in den Wasch-Port ( $\rightarrow$  Abb. 57, Seite 135) fährt:

**Hinweis:** Diese Prüfung sollten Sie vor der ersten Inbetriebnahme des Autosamplers durchführen.

1. Öffnen Sie den Riegel, mit dem Probennadel gesichert ist. Dadurch wird verhindert, dass sich die Probennadel verbiegt, wenn die Positionierung nicht korrekt sein sollte.



Abb. 48: Riegel am Nadelantrieb

2. Führen Sie einen Waschvorgang durch ( $\rightarrow$  Seite 120).

**Hinweis:** Vergewissern Sie sich, dass die Waschflasche mit Reinigungslösung gefüllt ist.

 Beobachten Sie, ob das Nadelpaar am Ende des Waschvorgangs korrekt in den Wasch-Port f\u00e4hrt. Dies ist der Fall, wenn sich die Nadel wie in Abb. 49 gezeigt im Waschport befindet und der Nadelhalter in der in Abb. 48 gezeigten Position verbleibt.



Abb. 49: Fused Silica-Nadel im Waschport

Ist der Nadelhalter erhöht, befindet sich das Nadelpaar nicht korrekt im Wasch-Port und muss neu justiert werden. Wenden Sie sich in diesem Fall bitte an den Kundendienst.

4. Stellen Sie sicher, dass die Nadel wieder richtig befestigt ist.

## 7.5.2 Waschen der Probennadel

Beim Waschen wird das Äußere der Probennadel im Waschport mit Waschflüssigkeit gespült und nach dem Waschen über die Membranpumpe mit Luft getrocknet. Es gibt zwei Möglichkeiten, die Probennadel zu waschen:

#### Über Chromeleon

- 1. Öffnen Sie in Chromeleon das Dialogfenster Commands für den Autosampler.
- 2. Legen Sie unter **WashVolume** das Volumen fest, das für den Waschvorgang aufgezogen werden soll.
- 3. Legen Sie unter **WashSpeed** die Geschwindigkeit der Spritze für den Waschvorgang fest.
- Legen Sie fest, über welche Position des Injektionsventils der Waschvorgang durchgeführt werden soll. Wählen Sie InjectValveToInject oder InjectValveToLoad. Bei der Einstellung InjectValveToLoad wird auch die Probenschleife mit gewaschen.
- 5. 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.

## 7.5.3 Wechseln der Probennadel

**Warnung:** Die Probennadel ist sehr spitz. Seien Sie vorsichtig, um sich nicht zu verletzen.

Beschreibung	BestNr.
WPS-3000(T)(B)PL (RS) Nano/Cap 2,4 µL-Probennadel, Fused Silica, nanoViper, schwarze Hülse	6820.3010
WPS-3000TBPL Analytisch und WPS-3000TBFC 15 µL-Probennadel, PEEK, Viper, rote Hülse	6820.3025
WPS-3000TFC 15 μL-Probennadel, Fused Silica, nanoViper, blaue Hülse	6820.3115
Informationen zu optional erhältlichen Probennadeln finden Sie im Kapitel 10.2 (→ Seite 171).	

1. Öffnen Sie den Riegel, mit dem die Probennadel gesichert ist.



Abb. 50: Riegel am Nadelantrieb

- 2. Geben Sie in Chromeleon das Kommando **PuncturerDownNeedleDown**. Damit werden der Vorstecher und die Probennadel nach unten gefahren.
- 3. Schrauben Sie die Probennadel am Injektions- bzw. Fraktionierventil ab (Port 6).
- 4. Hängen Sie die Probennadel aus der Halterung aus und ziehen Sie sie nach oben aus dem Vorstecher heraus.



Abb. 51: Entfernen der Probennadel

- 5. Führen Sie die neue Probennadel vorsichtig in den Vorstecher ( $\rightarrow$  Abb. 51) ein.
- 6. Hängen Sie die Probennadel in die Halterung ein ( $\rightarrow$  Abb. 51).
- 7. Schließen Sie die Probennadel an das Injektionsventil an ( $\rightarrow$  Abb. 7, Seite 28).
- 8. Schließen Sie den Riegel am Nadelantrieb ( $\rightarrow$  Abb. 50).
- 9. Führen Sie in Chromeleon das Kommando **PuncturerUpNeedleUp**. Damit werden der Vorstecher und die Probennadel nach oben bewegt.
- 10. Geben Sie das Kommando **SelfTest** über Chromeleon —*oder* rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Selftest**.
- Hinweis: Wenn die neue Nadel ein anderes Volumen hat als die Nadel, die Sie ausgebaut haben, sollten Sie sich vergewissern, dass in der Konfiguration des Autosamplers auf der Seite **Options** die Nadelgröße korrekt angegeben ist (→ Seite 44). Ändern Sie die Einstellung gegebenenfalls und übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Dazu steht Ihnen auf der Seite **General** die Schaltfläche **Send configuration to module** zur Verfügung (→ Seite 41).

## 7.5.4 Wechseln des Vorstechers

Warnung: Der Vorstecher ist sehr spitz. Seien Sie vorsichtig, um sich nicht zu verletzen.

Beschreibung	BestNr.
Vorstecher	6820.2400

- 1. Entfernen Sie die Probennadel wie in Kapitel 7.5.3 ( $\rightarrow$  Seite 121) beschrieben.
- Lösen Sie die Befestigungsschraube des Vorstechers mit dem Gabelschlüssel (Best.-Nr. 6000.0051) aus dem Gerätezubehör oder einem geschlitzten Ringschlüssel (Größe 1/4").



Lösen Sie die Befestigungsschraube für den Vorstecher.

Abb. 52: Befestigungsschraube Vorstecher

- 3. Führen Sie in Chromeleon das Kommando **PuncturerUpNeedleUp**. Damit werden der Vorstecher und die Probennadel nach oben bewegt.
- 4. Nehmen Sie die Schraube und den Vorstecher nach oben heraus.



Abb. 53: Vorstecher herausziehen

- 5. Setzen Sie einen neuen Vorstecher ein.
- 6. Schrauben Sie die Befestigungsschraube mit dem Sechskantschlüssel aus dem Gerätezubehör fest.
- 7. Installieren Sie die Probennadel wie in Kapitel 7.5.3 beschrieben.
- 8. Geben Sie das Kommando **SelfTest** über Chromeleon —*oder* rufen Sie am Gerät das Menü **Diagnostics** auf und wählen Sie **Selftest**.

Beschreibung	BestNr.
WPS-3000 (T)PL (RS) Nano/Cap und WPS-3000TBPL Nano/Cap 1 µL-Probenschleife, nanoViper, Fused Silica/PEEK sheathed	6826.2401
WPS-3000TFC 50 μL-Probenschleife, SST	6824.0019
WPS-3000TBPL Analytisch und WPS-3000TBFC 50 µL-Probenschleife, PEEK	6823.0019
Informationen zu optional erhältlichen Probenschleifen finden Sie im Kapitel 10.2 (→ Seite 171).	

## 7.6 Wechseln der Probenschleife

- 1. Schrauben Sie die Probenschleife am Injektionsventil ab.
- 2. Schalten Sie das Injektionsventil in die Position Load. Geben Sie dazu das Kommando InjectValveToLoad über Chromeleon —*oder* rufen Sie am Gerät das Menü Maintenance auf und wählen Sie Inject Valve.
- Führen Sie einen Waschvorgang mit mindestens 200 µL-Waschflüssigkeit bzw. mindestens dem doppelten Volumen der Probenschleife durch (→ Seite 120). Schließen Sie die neue Probenschleife an, während der Waschvorgang ausgeführt wird. Dies verhindert, dass Luft in das System gelangt.
- 4. Schließen Sie die Probenschleife am Port 2 des Injektionsventils an, sobald dort die Waschflüssigkeit austritt.
- 5. Schließen Sie die Probenschleife am Port 5 des Injektionsventils an, sobald die Waschflüssigkeit aus der Probenschleife austritt.
- **Hinweise:** Wenn die neue Probenschleife ein anderes Volumen hat als die Probenschleife, die Sie ausgebaut haben, sollten Sie sich vergewissern, dass
  - in der Konfiguration des Autosamplers auf der Seite Options

     (→ Seite 44) das Volumen der Probenschleife korrekt angegeben ist.
     Ändern Sie die Einstellung gegebenenfalls und übertragen Sie die neue Konfiguration von Chromeleon zum Autosampler. Dazu steht Ihnen auf der Seite General die Schaltfläche Send configuration to module zur Verfügung (→ Seite 41).
  - ♦ dass auch eine geeignete Pufferschleife und Spritze installiert sind (→ Seite 16).

## 7.7 Injektionsventil und Fraktionierventil

- **Vorsicht:** Um Schäden am Ventil zu vermeiden, achten Sie beim Betrieb des 900 bar-Ventils (WPS-3000(T)PL RS) darauf, das Ventil nicht trocken zu schalten.
- **Hinweis:** Um eine Verunreinigung der Ventilflächen zu vermeiden, tragen Sie Einmal-Handschuhe, wenn Sie Wartungsarbeiten am Ventil durchführen.

An zwei Ventilteilen, dem Rotor und dem Stator, kann es applikationsabhängig zu Verschleiß und damit zu Undichtigkeiten kommen. Beide Teile können unabhängig voneinander getauscht werden.

Ventile (2 Positionen, 6 Ports, 1/16")	BestNr.
WPS-3000(T)PL RS	
Rotor, SST (900 bar) Stator, SST (900 bar)	6826.0013 6826.0012
WPS-3000(T)PL Nano/Cap	
Rotor, SST Stator, SST	6820.0014 6820.0012
WPS-3000TBPL Nano/Cap und Analytisch	
Rotor, PAEK	6820.0035
Stator, PAEK	6820.0034
Stator, Titan	6820.0049
WPS-3000TFC	
Rotor Injektionsventil/Fraktionierventil, SST	6820.0014
Stator Injektionsventil/Fraktionierventil, SST	6820.0012
WPS-3000TBFC	
Rotor Injektionsventil/Fraktionierventil, PAEK	6820.0035
Stator Injektionsventil/Fraktionierventil, PAEK	6820.0034
Stator Injektionsventil, Titan	6820.0049

Fraktionierventil (2 Positionen, 6 Ports, 1/32")*	BestNr.
WPS-3000TBFC	
Rotor, PAEK Stator, PAEK	6720.0109 6720.0108

\*Nur zur Verwendung in einem nanoLC-System mit Tandem-Betrieb.

Fraktionierventile (2 Positionen, 8 Ports, 1/16")	BestNr.
WPS-3000(T)PL (RS) mit Micro Fraction Collection Option Rotor, SST Stator, SST	6820.0054 6820.0053
WPS-3000(T)BPL mit Micro Fraction Collection Option Rotor, PAEK Stator, PAEK	6821.0054 6821.0053

## 7.7.1 Reinigen des Ventils

Der erforderliche 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. Verwenden Sie eines der üblichen Lösungsmittel wie Methanol oder Acetonitril oder ein Gemisch (80/20) aus Methanol oder Acetonitril und Wasser.

## 7.7.2 Auseinanderbauen des Ventils

Auseinanderbauen eines 1/16"-Ventils



Abb. 54: Einzelteile des 1/16"-Ventils

- 1. Entfernen Sie mit einem Innensechskantschlüssel (Größe 9/64") die Statorschrauben. Lösen Sie abwechselnd jede Schraube jeweils nur um eine halbe Umdrehung, um ein Verkanten des Stators zu vermeiden.
- Entfernen Sie den Stator vom Ventilkörper. Legen Sie den Stator auf seine Außenseite, um die Dichtungsfläche nicht zu beschädigen. Sie können den Stator auch an den Kapillaren hängen lassen.
- 3. Entfernen Sie den Rotor vorsichtig per Hand oder mit einem kleinen Schraubendreher.
- 4. Überprüfen Sie die Dichtungsflächen des Rotors und Stators auf mögliche Kratzer:
  - Der Rotor bzw. Stator muss ausgetauscht werden, wenn Kratzer mit bloßem Auge erkennbar sind.
  - Wenn keine Kratzer erkennbar sind, reinigen Sie alle Teile gründlich mit einem geeigneten Lösungsmittel. Achten Sie darauf, die Oberfläche dabei nicht zu verkratzen. (Die häufigste Ursache für Probleme beim Einsatz des Ventils in der HPLC ist die Bildung von wasserlöslichen Pufferkristallen.) Nach der Reinigung muss der Rotor nicht getrocknet werden.

#### Auseinanderbauen eines 1/32"-Ventils

1. Entfernen Sie mit einem Innensechskantschlüssel (Größe 9/64") die Schrauben, die das Cap und den Stator am Ventil befestigen. Lösen Sie abwechselnd jede Schraube jeweils nur um eine halbe Umdrehung, um ein Verkanten des Stators zu vermeiden.



Abb. 55: Einzelteile des 1/32"-Ventils

- 2. Entfernen Sie die Ventilkappe. Eine polierte Dichtfläche befindet sich unten an der Ventilkappe. Am Stator befinden sich die Dichtflächen oben und unten. Um sicherzustellen, dass die Dichtflächen nicht beschädigt werden, legen Sie die Ventilkappe auf die Vorderseite der Kappe.
- 3. Entfernen Sie den Stator und legen Sie ihn auf eine saubere, weiche Oberfläche.
- 4. Entfernen Sie den Rotor vorsichtig per Hand oder einem kleinen Schraubendreher.
- 5. Prüfen Sie die Dichtflächen des Rotors und des Stators auf Kratzer.
  - Der Rotor bzw. Stator muss ausgetauscht werden, wenn Kratzer mit bloßem Auge erkennbar sind.
  - Wenn keine Kratzer erkennbar sind, reinigen Sie alle Teile gründlich mit einem geeigneten Lösungsmittel. Achten Sie darauf, die Oberfläche dabei nicht zu verkratzen. (Die häufigste Ursache für Probleme beim Einsatz des Ventils in der HPLC ist die Bildung von wasserlöslichen Pufferkristallen.) Nach der Reinigung muss der Rotor nicht getrocknet werden.

## 7.7.3 Zusammenbauen des Ventils

**i Hinweis:** Um eine Verunreinigung der Oberflächen zu vermeiden, reinigen Sie alle Ventilteile mit Isopropanol und tragen Sie Einmal-Handschuhe, wenn Sie das Ventil zusammenbauen.

#### Zusammenbauen eines 1/16"-Ventils

- 1. Setzen Sie den Rotor ein. Achten Sie auf die korrekte Ausrichtung: Die Dichtungsfläche muss mit den Flusswegen nach außen zeigen. Um zu verhindern, dass der Rotor versehentlich falsch eingesetzt wird, ist das Muster asymmetrisch.
- 2. Setzen Sie den Stator wieder auf. Ziehen Sie die beiden Schrauben handfest an. Ziehen Sie die Schrauben nicht zu fest an. Sie dienen lediglich der Befestigung und haben keinen Einfluss auf die Dichtkraft. Die Abdichtung erfolgt automatisch, wenn der Stator auf den Ventilkörper trifft.
- 3. Testen Sie das Ventil auf Dichtigkeit, indem Sie im System Druck herstellen. Wenn das Ventil nicht dicht ist, sollte es zur Reparatur eingeschickt werden.
- 4. Nach einem Rotor-Tausch

Führen Sie nach dem Austausch des Rotors in Chromeleon den Befehl **RotorSeal Changed** aus. Dadurch werden die entsprechenden Serviceinformationen aktualisiert.

 Folgen Sie den nachfolgenden Hinweisen nach dem Zusammenbau des Ventils im (→ nächste Seite).

#### Zusammenbauen eines 1/32"-Ventils

- 1. Setzen Sie den Rotor ein. Achten Sie auf die korrekte Ausrichtung: Die Dichtungsfläche muss mit den Flusswegen nach außen zeigen. Um zu verhindern, dass der Rotor versehentlich falsch eingesetzt wird, ist das Muster asymmetrisch.
- Setzen Sie den Stator wieder auf. Achten Sie auf die korrekte Ausrichtung: die Vorderseite muss vom Ventil weg zeigen. Vorder- und Rückseite des Stators unterscheiden sich darin, dass die Anschlüsse für die Kapillaren auf der Vorderseite größer sind und konisch kleiner werden bis zur Rückseite des Stators.
- 3. Setzen Sie die Ventilkappe wieder auf. Ziehen Sie die beiden Schrauben handfest an. Ziehen Sie die Schrauben nicht zu fest an. Sie dienen lediglich der Befestigung und haben keinen Einfluss auf die Dichtkraft. Die Abdichtung erfolgt automatisch, wenn der Stator auf den Ventilkörper trifft.
- 4. Testen Sie das Ventil auf Dichtigkeit, indem Sie im System Druck herstellen. Wenn das Ventil nicht dicht ist, sollte es zur Reparatur eingeschickt werden.

5. Nach einem Rotor-Tausch

Führen Sie nach dem Austausch des Rotors in Chromeleon den Befehl **RotorSeal Changed** aus. Dadurch werden die entsprechenden Serviceinformationen aktualisiert.

6. Folgen Sie den nachfolgenden Hinweisen nach dem Zusammenbau des Ventils.

#### Nach dem Zusammenbau des Ventils

Beachten Sie folgende Hinweise für den Anschluss von Kapillaren am Ventil:

- Wenn Sie mit einem biokompatiblen Autosampler arbeiten, beachten Sie bitte die Hinweise in Kapitel 8.3 (→ Seite 146).
- Verwenden Sie zum Anschluss der Kapillaren an das Ventil die vom Ventilhersteller empfohlenen Schneidringe und Fittingschrauben. Beachten Sie die Installationshinweise des Ventilherstellers.

## 7.8 Wechseln der Sicherungen



ng: Schalten Sie den Autosampler aus. Ziehen Sie den Netzstecker, bevor Sie die Sicherungen wechseln.

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



Abb. 56: Sicherungshalteschlitten

2. Tauschen Sie die Sicherungen aus.



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

Beschreibung	BestNr.
Autosampler <i>ohne</i> Probenthermostatisierung:	enthalten in: Kit WPS-Sicherungen
Sicherung, 2 A, träge, 5 x 20mm	(BestNr. 6820.0026)
Autosampler <i>mit</i> Probenthermostatisierung:	Informationen zum Inhalt des Kits finden Sie
Sicherung, 4 A, träge, 5 x 20mm	im Kapitel 10.3 (→ Seite 180).

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

## 7.9 Aktualisieren der Autosamplerfirmware

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

Welche Firmware-Version im Autosampler installiert und welche in Chromeleon enthalten ist, können Sie wie folgt feststellen:

- Firmware-Version des Autosamplers
  - Schalten Sie den Autosampler über den Netzschalter auf der Geräterückseite sein. Auf dem Autosamplerdisplay erscheinen allgemeine Informationen zum Autosampler, einschließlich der Firmware-Version.
  - Rufen Sie am Autosamplerdisplay das Menü **Diagnostics** auf (→ Seite 81) 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 dem nächsten Service Pack zu Chromeleon ausgeliefert und in den entsprechenden Release Notes beschrieben.

Die neue Firmware wird *nicht* automatisch auf den Autosampler übertragen, wenn Sie das Service Pack installieren. 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 zunächst, 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 40).
- 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 41) 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 bitte an den Kundendienst.

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

# 8 Autosampler-Spezifische Informationen

## 8.1 Funktionsprinzip WPS-3000PL (RS)

Die nachfolgende Abbildung illustriert das Funktionsprinzip des WPS-3000PL (RS):



Abb. 57: Funktionsprinzip

Der Autosampler arbeitet nach dem Pulled-Loop-Injektionsprinzip: Wenn das Injektionsventil noch in der Position **Inject** steht, wird die Probe über die Spritze aus dem Probenbehälter in die Pufferschleife gesaugt, so dass die Nadel vollständig mit Probe gefüllt ist. Danach wird das Injektionsventil in der Position **Load** umgeschaltet. Die Probe wird über die Probennadel in die Probenschleife gesaugt. Anschließend wird das Injektionsventil von der Position **Load** in die Position **Inject** umgeschaltet. Damit wird die Probe aus der Probenschleife über den Hochdruckflussweg auf die Trennsäule aufgebracht.

Die Probe wird im Back-Flush-Modus auf die Säule aufgebracht: In der Position Load wird die Probe in Richtung Spritze in die Probenschleife aufgezogen. In der Position Inject wird die Probe in entgegen gesetzter Richtung auf die Säule transportiert.



Abb. 58: Back-Flush-Modus

Der Autosampler enthält die folgenden Komponenten:

Komponente	Beschreibung
Injektionsventil / Probenschleife	Über das Injektionsventil und die Probenschleife wird die Probe in die mobile Phase eingebracht.
Nadelpaar	Das Nadelpaar besteht aus dem Vorstecher und der Probennadel. Der Vorstecher durchsticht vor der Injektion die Kappe/das Septum des Probenfläschchens bzw. die Versiegelung der Mikrotiterplatte. Über die Probennadel wird die Probe aufgezogen.
Wasch-Port	Im Wasch-Port können das Innere und Äußere der Probennadel gespült werden.
Drehteller	Der Drehteller ist mit den Mikrotiterplatten und/oder Probenhaltern bestückt.
Spritze	Über die Spritze wird die Probe aus der Kavität bzw. dem Probenfläschchen in die Probenschleife gesaugt.
Pufferschleife	Um die Spritze vor Verunreinigungen durch die Probe zu schützen, ist zwischen dem Injektionsventil und der Spritze eine Pufferschleife installiert.
Waschflüssigkeit	Mit Hilfe der Waschflüssigkeit wird die restliche Probe aus der Pufferschleife und der Probennadel entfernt sowie die Schleife und die Probennadel gespült.
Luftpumpe	Mit Hilfe der Luftpumpe kann die Außenseite der Probennadel getrocknet werden.
X, Z und z'	Bezeichnen die möglichen Bewegungsrichtungen des Nadelpaars.
Y	Bezeichnet die mögliche Bewegungsrichtung des Drehtellers.

Informationen zu den unterstützten Injektionsmethoden finden Sie im technischen Anhang ( $\rightarrow$  Seite 150).

Nahezu alle Funktionen stehen über benutzerdefinierte Programme (UDP = user-defined program) in Chromeleon zur Verfügung ( $\rightarrow$  Seite 96).
# 8.2 Autosampler zur Fraktionssammlung

# 8.2.1 WPS-3000FC

Der WPS-3000FC ist mit *zwei* Ventilen ausgestattet, einem Injektionsventil und einem Fraktionierventil, so dass Injektion, Fraktionierung und erneute Injektion vollautomatisch auf einem einzigen System durchgeführt werden können.



Abb. 59: WPS-3000TFC (hier biokompatible Konfiguration)

Das Funktionsprinzip des Autosamplers ist in Kapitel 8.2.1.1 (→ Seite 138) kurz erläutert.

Den Autosampler gibt es auch in einer biokompatiblen Version. Beachten Sie dazu bitte die Hinweise zum Anschluss der Komponenten an den Ventilen ( $\rightarrow$  Seite 146). Das biokompatible Fraktionierventil kann durch ein Ventil mit 0,1 mm Bohrung (1/32") ersetzt werden (Best.-Nr. 6825.0020).

Das Injektionsventil kann gegen ein 900 bar-Ventil ausgetauscht werden. Wenden Sie sich dazu bitte an den Kundendienst.

# Umrüstkits

#### Nano/Kap-Umrüstkit für Nano/Kap-Anwendungen

Zusätzlich steht ein Umrüstkit zur Verfügung, mit denen der Autosampler für Anwendungen in der Nano/Kap-LC optimiert werden kann ( $\rightarrow$  Seite 143).

#### Normalphasen-Kit für WPS-3000TFC Autosampler

Ein Normalphasen Umrüstkit steht zur Umrüstung des WPS-3000TFC Autosamplers für Normalphasen-Anwendungen zur Verfügung. Informationen zum Kit finden Sie in Kapitel 8.2.1.5 ( $\rightarrow$  Seite 144).

# 8.2.1.1 Funktionsprinzip

Die nachfolgende Abbildung illustriert das Funktionsprinzip des WPS-3000FC:



Abb. 60: Funktionsprinzip

Der Autosampler arbeitet nach dem Pulled-Loop-Injektionsprinzip: Wenn das Injektionsventil in der Position **Inject** und das Fraktionierventil in der Position **Drain** steht, wird die Probe über die Spritze aus dem Probenbehälter in die Pufferschleife gesaugt, so dass die Nadel und die Verbindungskapillare zwischen den beiden Ventilen (Bridge) vollständig mit Probe gefüllt ist. Danach wird das Injektionsventil in der Position **Load** umgeschaltet. Die Probe wird über die Probennadel und die Bridge in die Probenschleife gesaugt. Anschließend wird das Injektionsventil von der Position **Load** in die Position **Inject** umgeschaltet. Damit wird die Probe aus der Probenschleife über den Hochdruckflussweg auf die Trennsäule aufgebracht.

Die Probe wird im Back-Flush-Modus auf die Säule aufgebracht: In der Position Load wird die Probe in Richtung Spritze in die Probenschleife aufgezogen. In der Position Inject wird die Probe in entgegen gesetzter Richtung auf die Säule transportiert.

Nach der Injektion kann das Fraktionierventil in die Position **Collect** geschaltet werden, um über die Probennadel Fraktionen in die Probenbehälter zu sammeln.

Komponente	Beschreibung
Injektionsventil / Probenschleife	Über das Injektionsventil und die Probenschleife wird die Probe in die mobile Phase eingebracht.
Fraktionierventil	Wenn das Fraktionierventil nach der Injektion in die Position <b>Collect</b> geschaltet wird, können über die Probennadel Fraktionen in die Probenbehälter gesammelt werden.
Bridge	Bezeichnet die Verbindungskapillare zwischen Injektions- und Fraktionierventil.
Nadelpaar	Das Nadelpaar besteht aus dem Vorstecher und der Probennadel. Der Vorstecher durchsticht vor der Injektion die Kappe/das Septum des Probenfläschchens bzw. die Versiegelung der Mikrotiterplatte. Über die Probennadel wird die Probe aufgezogen.
Wasch-Port	Im Wasch-Port können das Innere und Äußere der Probennadel gespült werden.
Drehteller	Der Drehteller ist mit den Mikrotiterplatten und/oder Probenhaltern bestückt.
Spritze	Über die Spritze wird die Probe aus der Kavität bzw. dem Probenfläschchen in die Probenschleife gesaugt.
Pufferschleife	Um die Spritze vor Verunreinigungen durch die Probe zu schützen, ist zwischen dem Injektionsventil und der Spritze eine Pufferschleife installiert.
Waschflüssigkeit	Mit Hilfe der Waschflüssigkeit wird die restliche Probe aus der Pufferschleife und der Probennadel entfernt sowie die Schleife und die Probennadel gespült.
Luftpumpe	Mit Hilfe der Luftpumpe kann die Außenseite der Probennadel getrocknet werden.
X, Z und z'	Bezeichnen die möglichen Bewegungsrichtungen des Nadelpaars.
Y	Bezeichnet die mögliche Bewegungsrichtung des Drehtellers.

Der Autosampler enthält die folgenden Komponenten:

Informationen zu den unterstützten Injektionsmethoden finden Sie im technischen Anhang ( $\rightarrow$  Seite 150).

Nahezu alle Funktionen stehen über benutzerdefinierte Programme (UDP = user-defined program) in Chromeleon zur Verfügung ( $\rightarrow$  Seite 96).

#### 8.2.1.2 Injektions- und Fraktionierventile

Der Autosampler ist mit einem Injektionsventil und einem Fraktionierventil (je 2 Positionen, 6 Ports) ausgestattet. Das Fraktionierventil ist geeignet für Anwendungen mit folgendem Druck (Angaben zum Injektionsventil siehe Seite 28):

Fraktionierventil	Material	Maximaler Druck	Beschreibung
WPS-3000TFC	Edelstahl	3,5 MPa (35 bar)	1/16", 0,25 mm, 75°C
WPS-3000TBFC	PAEK	3,5 MPa (35 bar)	1/16", 0,25 mm, 50°C
WPS-3000TBFC (optional bestellbar)	PAEK	35 MPa (350 bar)	1/32", 0,1 mm, 50°C

Jedes Ventil hat zwei Betriebspositionen. Das Injektionsventil hat die Positionen Load und Inject, das Fraktionierventil die Positionen Collect und Drain.

Für eine Injektion müssen das Injektionsventil in der Position **Load** und das Fraktionierventil in der Position **Drain** stehen. Die Probe dann wird über die Probennadel und Verbindungskapillare zwischen den beiden Ventilen (Bridge) in die Probenschleife gesaugt. Wenn das Injektionsventil in die Position **Inject** umgeschaltet, wird die Probe in den Hochdruckflussweg zur Analyse auf die Trennsäule aufgebracht.

Nach der Injektion kann das Fraktionierventil in die Position **Collect** geschaltet werden, um Fraktionen über die Probennadel in die Probenfläschchen zu sammeln.



Die Komponenten sind wie folgt an den Ventilen angeschlossen:

Abb. 61: Anschlüsse am Injektions- und Fraktionierventil

Injektionsventil			
Port Anschluss von			
1	Spritze/Pufferschleife		
2	Probenschleife		
3	Pumpe		
4	Säule		
5	Probenschleife		
6	Verbindungskapillare (Bridge)		

F Port	raktionierventil Anschluss von
1	Verbindungskapillare (Bridge)
2	Abfallleitung
3	Abfallleitung
4	Abfall
5	Säule/Detektor
6	Probennadel

Wenn Sie mit einem biokompatiblen Autosampler arbeiten, beachten Sie bitte auch die Hinweise zum Anschluss der Kapillaren an den Ventilen ( $\rightarrow$  Seite 146).

# 8.2.1.3 Hinweise zum Betrieb

Beachten Sie die folgenden Hinweise für den Betrieb des Autosamplers:

- Das Sammeln von Fraktionen wird in Chromeleon 7.1 oder später derzeit noch nicht unterstützt.
- Zur Steuerung des Autosamplers unter Chromeleon werden Chromeleon 6.80 Service Release 6 (oder höher) und folgende Lizenzen benötigt:
  - Timebase Class 1
  - Fraction Collection (Best.-Nr. 5960.0038) mit den *Grundfunktionen* f
    ür die Fraktionssammlung, mit Programmassistent und Fraktionsreport.

oder

• Extended Fraction Collection (Best.-Nr. 5960.0039)

für die *erweiterte* Fraktionssammlung. Die Lizenz **Extended Fraction Collection** schließt die Lizenz **Fraction Collection** ein und unterstützt zusätzlich fortgeschrittene Funktionen zur Fraktionssammlung wie Automatische Aufreinigung von Proben über die zugehörigen Post-acquisition steps, Zuordnung von Proben und Fraktionen anhand von Farben, ausgefeilten Fraktionieralgorithmen mit Peakschulter-Erkennung

Weitere Informationen zu den beiden Lizenzen erhalten Sie von der an den Thermo Fisher Scientific Vertriebsorganisation für Dionex HPLC-Produkte.

 Vergewissern Sie sich, dass in der Konfiguration des Autosamplers auf der Seite Options (→ Seite 45) das Kontrollkästchen WPS-3000FC/WPS-3000TFC aktiviert und das BridgeVolume korrekt angegeben ist.

Das **BridgeVolume** ist das Volumen der Kapillare, die das Injektionsventil mit dem Fraktionierventil verbindet. Vergewissern Sie sich dass, der angegebenen Wert mit dem Wert auf dem Label der Kapillare übereinstimmt. Ist das BridgeVolume nicht korrekt angegeben, kann es zu Fehlern bei der Injektionssteuerung kommen.

- **I** Hinweis: Denken Sie daran, den Wert im Feld **BridgeVolume** entsprechend anzupassen, wenn Sie zwischen den beiden Ventilen eine Kapillare mit einem anderen Volumen installieren oder wenn Sie den Autosampler für eine Applikation einsetzen, bei der die beiden Ventile *nicht* miteinander verbunden sind (in diesem Fall ist BridgeVolume auf 0 zu setzen).
- Damit die Funktionen zur Fraktionssammlung in Chromeleon verfügbar sind, muss zusätzlich zum Autosampler der Treiber **Fraction Collection** installiert werden:
  - a) Installieren Sie den Treiber in der Zeitbasis, in der auch der Autosampler installiert ist. Gehen Sie analog der Beschreibung zur Installation des Autosamplers vor (→ Seite 40).

b) Wählen Sie unter **Manufacturers** den Eintrag **Generic** und unter **Devices** den Eintrag **Fraction Collection** aus.

Auf der Seite **General** werden unter **Device Name** der Name, unter dem der Treiber in der Installationsumgebung und im Chromeleon Client geführt wird, und die maximale Anzahl an Detektionskanälen angezeigt. Prüfen und ändern Sie die Angaben, falls erforderlich.

Wenn Sie vorhandene Steuerfenster verwenden möchten, sollten Sie den vorgegebenen **Device Name** nicht ändern.

- c) Klicken Sie **Ok**, um die Installation zu abzuschließen.
- d) Speichern Sie die Konfiguration mit **Save Installation** im Menü **File** und schließen Sie das Konfigurationsprogramm.
- Ab Chromeleon 6.80 SR10 steht eine verbesserte Unterstützung der Fraktionierung durch den Programmassistenten und spezielle Steuerfenster zur Verfügung. Beim Erstellen eines Programms mit dem Programmassistenten werden alle Programmteile für das Fraktionieren erzeugt ein manuelles Nachbearbeiten ist nicht erforderlich. Dazu gibt es zwei zusätzliche Wizard-Seiten, auf denen Sie die erste Fraktionierposition, das Meanderschema, die Airdry-Option und die Optionen zu den Fraktionsgefäßen einstellen können.
  - Fraktionierprogramme sind spezifisch für einen bestimmten Fraktionenhalter. Wenn Sie einen anderen Fraktionenhalter konfigurieren, müssen Sie auch den Fraktionenhalter auf der Seite **Collector General Options** des PGM-Editors ändern.
  - Parameter, die das Tray-Layout und das Fraktionierschema sowie die Nummerierung der Fraktionen betreffen dürfen während des gesamten Batches nicht geändert werden.



*Abb.* 62: *Abschnitt zur Fraktionierung auf dem Steuerfenster des WPS-3000FC* 

Weitere Informationen zur Fraktionssammlung und eine Liste der für den Autosampler verfügbaren Kommandos und Properties finden Sie in der *Chromeleon Hilfe*.

# 8.2.1.4 Optimieren des Autosamplers für Nano/Cap-Anwendungen

Beide Autosamplerversionen können für Anwendungen in der Nano- und Cap-LC optimiert werden.

Beschreibung	BestNr.
Umrüstkit Nano/Kap für WPS-3000TFC mit - 2,4 μL-Probennadel (Fused Silica, nanoViper) - 25 μL-Spritze - 50 μL-Pufferschleife - 1 μL-Probenschleife, nanoViper - 1 μL-Bridge (PEEKsil <sup>TM</sup> )	6824.0030
<i>Optional erhältliche Probenschleifen (nanoViper)</i> : 5 μL 10 μL 20 μL	6826.2405 6826.2410 6826.2420
Umrüstkit Nano/Kap für WPS-3000TBFC mit - 2,4 μL-Probennadel (Fused Silica, nanoViper) - 25 μL-Spritze - 50 μL-Pufferschleife - 1 μL-Probenschleife, nanoViper - 1 μL-Bridge (PEEKsil, biokompatibel)	6825.0030
<i>Optional erhältliche Probenschleifen (nanoViper)</i> : 5 μL 10 μL 20 μL	6826.2405 6826.2410 6826.2420

Installieren Sie die Komponenten aus dem Umrüstkit wie in den jeweiligen Servicekapiteln beschrieben ( $\rightarrow$  Kapitel 7, ab Seite 111 und 8.2.1.6, Seite 144).

# 8.2.1.5 Optimieren des WPS-3000TFC für Normalphasen-Anwendungen

Ein Normalphasen-Kit (Best.-Nr. 6820.0060) steht **nur** für den WPS-3000TFC Autosampler (Best.-Nr. 5824.0020) zur Verfügung. Verwenden Sie Normalphasen (NP)-Eluenten **nur** mit dem WPS-3000TFC Autosampler, nachdem der Autosampler mit dem Normalphasen-Kit umgerüstet wurde.

# **STOP** Warnung

Beachten Sie für NP-Applikationen die Sicherheitshinweise zu den WPS-3000TFC und WPS-3000TBFC Autosamplern auf Seite 5.

Beschreibung	BestNr.
Normalphasen-Kit, enthält 1 Nadel, 15 µL, SST, Viper (200 µm x 0,5 mm (I.D. x A.D.)), mit Erdungskabel 1 Kapillare, Viper (SST, 0,18 x 250 mm (I.D. x L), zur Verwendung als 6,4 µL-Bridge) 1 Sachskent Stiffachläggel (Cräße 2 mm)	6820.0060
1 Unterlegscheibe und 1 Sicherungsscheibe	

# 8.2.1.6 Service

Die Servicemaßnahmen und Reparaturarbeiten für den Autosampler sind im Kapitel Service beschrieben ( $\rightarrow$  ab Seite 111).

Zusätzlich kann beim WPS-3000TFC bei Bedarf die Verbindungskapillare (Bridge) zwischen Injektionsventil und Fraktionierventil getauscht werden.

Beschreibung	BestNr.
Bridge (PEEK, Volumen: 6,2 μL) für WPS-3000TFC WPS-3000TBFC	6824.0016 6825.0016
<i>Bridge (PEEKsil, Volumen: 1 μL) für</i> WPS-3000TFC, optimiert für Nano/Kap-Anwendungen WPS-3000TBFC, optimiert für Nano/Kap-Anwendungen	6824.0015 6825.0015

Lösen Sie die Kapillare am Port 6 des Injektionsventils und am Port 1 des Fraktionierventils. Wenn Sie eine Kapillare durch eine Kapillare mit einem anderen Volumen ersetzen, müssen Sie den Wert für **BridgeVolume** in den Eigenschaften des Autosamplers entsprechend anpassen (→ Seite 45). Andernfalls kann es zu Fehlern bei der Injektionssteuerung kommen.

# 8.2.2 Autosampler mit Micro Fraction Collection (MFC) Option

Für die Nano- und Kap-Autosampler WPS-3000(T)PL und WPS-3000TBPL bietet an den Thermo Fisher Scientific eine Micro Fraction Collection (MFC)-Option zum Sammeln von Fraktionen an. Mit dieser Option können Injektion, Fraktionierung und erneute Injektion vollautomatisch auf einem einzigen System durchgeführt werden.

Der Autosampler ist mit *einem* Injektions- und Fraktionierventil (2 Positionen, 8 Ports) ausgestattet. Das Ventil ist geeignet für Anwendungen mit folgendem Druck:

Ventil	Material	Maximaler Druck	Beschreibung
WPS-3000PL mit MFC	Edelstahl	35 MPa (350 bar)	1/16", 0,25 mm, 75°C
WPS-3000BPL mit MFC	PAEK	35 MPa (350 bar)	1/16", 0,25 mm, 50°C

Vorsicht: WPS-3000(T)PL und WPS-3000TBPL Nano/Cap-Autosampler mit MFC dürfen nicht für Normalphasen-Anwendungen eingesetzt werden.

Detaillierte Informationen zu dieser Option finden Sie in den begleitenden Dokumenten *Micro-Fraction Collection Option for the UltiMate 3000 Series Nano/Cap Autosampler* (in englischer Sprache):

- Installation Instructions mit Informationen zur Installation und Qualifizierung
- *Operating Instructions* mit Informationen zu Programmdateien in Chromeleon und Beispielen

# 8.3 Biokompatible Autosampler

Dieses Kapitel bezieht sich auf biokompatible Autosampler, bei denen das Injektionsventil (bzw. das Injektionsventil und Fraktionierventil) in PAEK ausgeführt ist.

Die am Ventil angeschlossenen Komponenten (Pufferschleife, Probennadel, Probenschleife, Säule, Kapillaren) bestehen aus PEEKsil, PEEK, PTFE oder Fused Silica. Da die Anschlussports bei PAEK-Ventilen tiefer ausgeführt sind ( $\rightarrow$  Abb. 63), müssen die Fittinge bei herkömmlichen Fittingverbindungen wie unten beschrieben bis zum Anschlag in den Port montiert sein, damit ein optimaler Anschluss ohne Totvolumen gewährleistet ist. Dies ist nicht notwendig für nanoViper- bzw. Viper-Kapillarverbindungen, die als totvolumenfreie Verbindungen konzipiert sind.



Abb. 63: Anschluss der Kapillaren am Injektionsventil

Verwenden Sie zum Anschluss das Werkzeug zur Fittingvormontage ( $\rightarrow$  Abb. 65), um Schäden am Ventil oder den Komponenten zu vermeiden. Das Werkzeug ist im Zubehör (WPS-3000TBPL Analytisch und WPS-3000TBFC) oder im Applikationskit (WPS-3000TBPL) enthalten und ist auch separat erhältlich (Best.-Nr. 6000.0065).

▲ Vorsicht: Verwenden Sie zum Anschluss der Komponenten am PAEK-Ventil keine Verschraubungen und/oder Schneidringe aus Edelstahl. Um eine Beschädigung des Ventils zu vermeiden, sollten Sie ausschließlich die mitgelieferten Kapillaren und Fittinge sowie original Dionex-Ersatzteile verwenden.

Schließen Sie die Komponenten wie unten beschrieben am biokompatiblen Schaltventil an. Die folgende Beschreibung beschreibt den Anschluss einer Nano-Säule, gilt aber analog für die anderen Komponenten. 1. Für eine totvolumenfreie Fittingverbindung zwischen der Säule und dem Schaltventil werden eine fingerfeste Verschraubung (finger-tight nut), ein Schneidring (ferrule) und ein Sleeve benötigt. Schieben Sie die Verschraubung und den Schneidring wie in Abb. 64 gezeigt auf den Sleeve.



Abb. 64: Fittingverbindung

2. Schieben Sie den Sleeve mit dem Schneidring und der Verschraubung in einen der Ports des Werkzeugs für die Fittingmontage und ziehen Sie die Verschraubung 2 bis 3 Umdrehungen mit der Hand an.



Abb. 65: Werkzeug zur Fittingvormontage

- ▲ Vorsicht: Um Schäden am biokompatiblen Schaltventil zu vermeiden, darf die Fittingverbindung keinesfalls am Schaltventil vormontiert werden. Ebenso sollte zum Anziehen der fingerfesten Verschraubung kein Werkzeug verwendet werden.
- 3. Schieben Sie den Sleeve bis zum Anschlag in den Port. Dies ist für eine totvolumenfreie Verbindung unerlässlich.



Abb. 66: Vormontage der Fittingverbindung

4. Drehen Sie die Verschraubung handfest in den Port hinein.

- 5. Ziehen Sie die Verschraubung vorsichtig eine weitere Viertelumdrehung (90°) über den Punkt hinaus fest, an dem der Schneidring den Sleeve zu fassen beginnt.
  - **Hinweis:** Eine Angabe des Drehmoments hinsichtlich der zum Anziehen der Verschraubung erforderlichen Kraft ist an dieser Stelle nicht möglich. Die Kraft kann aufgrund der Reibung zwischen der Verschraubung und dem Gewinde sowie des Materials und der Wandstärke der anzuschließenden Komponente variieren.
- 6. Nehmen Sie das vormontierte Fitting aus dem Port heraus und überprüfen Sie es. Der Schneidring sollte fest auf dem Sleeve installiert sein und sich nicht bewegen, wenn Sie vorsichtig daran ziehen. Wenn der Schneidring noch seitlich bewegt werden kann (→ Abb. 67), schieben Sie das Fitting nochmals in das Werkzeug hinein und ziehen Sie die Verschraubung eine weitere 1/8-Umdrehung fest (über handfest hinaus).



Abb. 67: Prüfen der Fittingverbindung

- 7. Entnehmen und prüfen Sie das Fitting. Wiederholen Sie den vorherigen Schritt falls erforderlich.
- 8. Schieben Sie die Säule (bzw. die Fused-Silica-Kapillare) bis zum Anschlag durch das Sleeve, bis diese aus dem Sleeve herausragt.



Abb. 68: Einführen der Säule

9. Schieben Sie die Säule zurück und befestigen Sie das vormontierte Fitting mit der Säule am entsprechenden Port des Schaltventils. Stellen Sie sicher, dass die Säule beim Befestigen am Port nicht aus dem Sleeve herausragt.



Abb. 69: Zurückschieben der Säule

- 10. Ziehen Sie die Verschraubung fingerfest an. Schieben Sie dabei Säule und Sleeve vorsichtig in das Ventil, so dass das Ende der Säule bis zum Anschlag in den Ventilport geschoben wird und kein Totvolumen entsteht.
- 11. Ziehen Sie die Verschraubung um eine weitere ¼-Umdrehung fest (über fingerfest hinaus).
- 12. Starten Sie den Fluss und prüfen Sie, ob die Verbindung dicht ist. Stoppen Sie bei Undichtigkeiten den Fluss und lösen Sie das Fitting. Wiederholen Sie die Schritte 8 bis 11.
- I Hinweis: Beim WPS-3000TBPL sind in den Applikationskits für das biokompatible UltiMate 3000-System lange Hex-Verschraubungen (→ Abb. 70; Best.-Nr. 6721.0017) enthalten, die Sie anstatt der beschriebenen Fittingverbindungen verwenden können. Schließen Sie die Hex-Verschraubung an den entsprechenden Port des Injektionsventils an und ziehen Sie die Verschraubung mit Hilfe eines Sechskantschlüssels (Größe: ¼") fest.



Abb. 70: Lange Hex-Verschraubung

# 8.4 Injektionsmethoden

Der Autosampler unterstützt die folgenden Injektionsmethoden. Informationen dazu, wie die Injektionsmethode in Chromeleon festgelegt werden kann, finden Sie auf Seite 83. Informationen zu *benutzerdefinierten* Injektionsprogrammen finden Sie auf Seite 96 und in der Hilfe zu Chromeleon.

Die Injektionsmethoden sind nachfolgend am Beispiel des WPS-3000PL beschrieben.

# 8.4.1 Full-Loop-Injektionen

Full-Loop-Injektionen bieten höchste Reproduzierbarkeit (RSD < 0,4% bei 1µL-Injektionen), sind aber nicht ganz so genau. Der Probenbedarf berechnet sich wie folgt:

Für die erste Injektion:

FlushVolume	+	LoopVolume	x	LoopOverfill	
Beispiel:					
5,0 μL	+	5,0 μL	Х	2	= 15 µL

Für alle weiteren Injektionen aus demselben Probengefäß:

FlushVolume2	+	LoopVolume	x	LoopOverfill	
Beispiel:					
2,4 μL	+	5,0 μL	х	2	= 12,4 μL

**Hinweis:** Bei Verwendung der Fused-Silica-Probennadel sollte das Spülvolumen mindestens 5.0 µL betragen. Bei hochviskosen Proben kann es unter Umständen erforderlich sein, zur Verbesserung der Genauigkeit das Spülvolumen zu erhöhen und eine langsamere Spritzengeschwindigkeit zu wählen.

Nachfolgend ist die Schaltsequenz für eine Standard-Full-Loop-Injektion beschrieben. Ergänzend dazu unterstützt der Autosampler eine Optimierung des Injektionsprofils ( $\rightarrow$  Seite 160).

Zu Beginn der Sequenz befindet sich das Ventil in der Position INJECT ( $\rightarrow$  Abb. 71). Der Vorstecher hat die Versiegelung/das Septum des Probengefäßes durchstoßen und die Probennadel befindet sich im Probengefäß.



Abb. 71: Position INJECT

Die Spritze zicht das programmierte FlushVolume aus dem Probengefäß auf. Indem die Probennadel mit Probe gefüllt wird ( $\rightarrow$  Abb. 72), wird die restliche Waschflüssigkeit, die sich eventuell noch in der Kapillare befindet, daraus entfernt.



Abb. 72: Aufziehen von Probe in die Probennadel

Wenn das Injektionsventil in die Position LOAD schaltet ( $\rightarrow$  Abb. 73), hat die Probe am Einlass in die Probenschleife eine "scharfe" Kante.



Abb. 73: Position LOAD

Bei Full-Loop-Injektionen wird die Probenschleife mengenmäßig befüllt, indem ein Vielfaches (definiert durch den LoopOverfill-Faktor) des Schleifenvolumens durch die Schleife transportiert wird ( $\rightarrow$  Abb. 74).



Abb. 74: Transport des Schleifenvolumens durch die Probenschleife

Das Injektionsventil schaltet in die Position INJECT ( $\rightarrow$  Abb. 75). Die Probenschleife wird Teil des Flussweges der mobilen Phase; die Probe wird auf die Säule transportiert und die Analysenzeit beginnt.



Abb. 75: Position INJECT

Die nachfolgende Abbildung zeigt die Verteilung der Flüssigkeiten bei einer Full-Loop-Injektion kurz vor der Injektion (wenn zuvor ein Wash ausgeführt wurde):

Nadel <sub>i</sub> Probenschleife	Pufferschleife				
					$\Box$
Probe		Eluent	Probe	Spülflüssigkeit	

Abb. 76: Full-Loop-Injektion

**i** Hinweise: Wenn aus einem Probengefäß nur ein Mal injiziert wird oder wenn nach jeder Injektion ein Waschzyklus durchgeführt werden muss, wird die Probennadel direkt nach der Injektion aus dem Probengefäß entfernt (und sofort gewaschen, sofern ein Waschzyklus programmiert wurde). Nach Ablauf der Analysenzeit wird eine neue Probe gestartet.

Wenn aus einem Probengefäß mehrfach injiziert wird, ohne dass ein Waschzyklus programmiert wurde, wird das programmierte FlushVolume2 verwendet. Wenn die Probenmenge, die für die nächste Injektion aus dem Probengefäß aufgezogen wird, das Gesamtvolumen der Pufferschleife übersteigt, wird vor der nächsten Injektion ein Waschvorgang (Wash) durchgeführt. Zu Beginn der nächsten Probe wird dann wieder das gesamte programmierte Spülvolumen (FlushVolume) aufgezogen.

Der Wert für FlushVolume2 kann kleiner sein als der für FlushVolume. Thermo Fisher Scientific empfiehlt, den Wert immer mindestens gleich oder größer zu wählen als das Nadelvolumen.

# 8.4.2 Partial-Loop-Injektionen

Partial-Loop-Injektionen bieten höchste Genauigkeit. (Bitte beachten Sie dabei, dass die Genauigkeit von der Genauigkeit der Spritze abhängt.) Bei Injektionsvolumen > 200 nL, ist die Reproduzierbarkeit besser als 1% RSD. Empfohlen wird ein Spülvolumen von 5,0  $\mu$ L. Sie können auch ein niedrigeres Spülvolumen wählen. Dies kann jedoch die Leistungsfähigkeit beeinträchtigen.

Im Hinblick auf beste Reproduzierbarkeit und Genauigkeit wird die Verwendung einer 25  $\mu$ L-Spritze mit einer 1 $\mu$ L-Probenschleife empfohlen. Damit vermeiden Sie einen Genauigkeitsverlust aufgrund der Expansion des Schleifeninhalts, wenn das Injektionsventil vor dem Aufziehen der Probe von der Position **INJECT** in die Position **LOAD** geschaltet wird. Bei einem hohen Druck (200 bar), kann dieser Verlust für eine 5  $\mu$ L-Schleife bis zu 0,025  $\mu$ L betragen.

Für Partial-Loop-Injektionen berechnet sich der erforderliche Mindest-Probenbedarf wie folgt:

Für die erste Injektion:

FlushVolume + Injektionsvolumen Beispiel:  $5,0 \ \mu L$  +  $2,5 \ \mu L$  =  $7,5 \ \mu L$ 

Für alle weiteren Injektionen aus demselben Probengefäß:

FlushVolume2	+	Injektionsvolumen	
Beispiel:			

 $2,4 \ \mu L$  +  $2,5 \ \mu L$  =  $4,9 \ \mu L$ 

Wenn zwischen den Injektionen ein Waschzyklus durchgeführt wird, ist der Probenverlust immer das FlushVolume. Wählen Sie für Injektionen ohne Probenverlust den Mikroliter-Pickup-Modus

Nachfolgend ist die Schaltsequenz für eine Standard-Partial-Loop-Injektion beschrieben. Ergänzend dazu unterstützt der Autosampler eine Optimierung des Injektionsprofils ( $\rightarrow$  Seite 160).

Die ersten drei Schritte entsprechen denen bei einer Full-Loop-Injektion ( $\rightarrow$  Seite 150).

Bei Partial-Loop-Injektionen wird das programmierte Injektionsvolumen in die Probenschleife transportiert ( $\rightarrow$  Abb. 77).



Abb. 77:Befüllen der Probenschleife mit dem gewünschten Injektionsvolumen

**I** Hinweis: Das Injektionsvolumen kann frei gewählt werden. Wenn das Injektionsvolumen 50% des Probenschleifenvolumens übersteigt, nimmt die Linearität stark ab. In Chromeleon wird dann eine Warnung ausgegeben. Wenn das Injektionsvolumen gleich dem Probenschleifenvolumen ist, wird ein Error ausgegeben.

Das Injektionsventil schaltet in die Position INJECT ( $\rightarrow$  Abb. 78). Die Probenschleife wird Teil des Flussweges der mobilen Phase; die Probe wird auf die Säule transportiert und die Analysenzeit beginnt.



Abb. 78: Position INJECT

Die folgende Abbildung zeigt die Verteilung der Flüssigkeiten bei einer Partial-Loop-Injektion:



Abb. 79: Partial-Loop-Injektion

Hinweise: Wenn aus einem Probengefäß nur ein Mal injiziert wird oder wenn nach jeder Injektion ein Waschzyklus durchgeführt werden muss, wird die Probennadel direkt nach der Injektion aus dem Probengefäß entfernt (und sofort gewaschen, sofern ein Waschzyklus programmiert wurde). Nach Ablauf der Analysenzeit wird eine neue Probe gestartet.

Wenn aus einem Probengefäß mehrfach injiziert wird, ohne dass ein Waschzyklus programmiert wurde, wird das programmierte FlushVolume2 verwendet. Wenn die Probenmenge, die für die nächste Injektion aus dem Probengefäß aufgezogen wird, das Gesamtvolumen der Pufferschleife übersteigt, wird vor der nächsten Injektion ein Waschvorgang (Wash) durchgeführt. Zu Beginn der nächsten Probe wird dann wieder das gesamte programmierte Spülvolumen (FlushVolume) aufgezogen.

Der Wert für FlushVolume2 kann kleiner sein als der für FlushVolume. Thermo Fisher Scientific empfiehlt, den Wert immer mindestens gleich oder größer zu wählen als das Nadelvolumen.

# 8.4.3 Mikroliter Pick-up

Mikroliter-Pick-up erzeugt keinen Probenverlust und höchste Genauigkeit, aber eine etwas geringere Reproduzierbarkeit. Verwenden Sie eine 25µL-Spritze und eine 10µL-Probenschleife, um eine optimale Genauigkeit und Reproduzierbarkeit zu erreichen. Es wird ein Pfropfen Transportflüssigkeit, gefolgt von dem Probenpfropfen in die Probenschleife transportiert, wiederum gefolgt von Transportflüssigkeit. Das Volumen des Transportflüssigkeitspfropfens kann über das Property FlushVolume festgelegt werden.

Um die Genauigkeit beim Mikroliter Pick-up zu gewährleisten, gibt es eine Obergrenze für die injizierbare Probenmenge. Diese Grenze, die sich aus den Volumen der Probenschleife und der Probennadel ergibt, kann wie folgt berechnet werden:

#### 1/3 LoopVolume

Bei Verwendung einer 20-µL Probenschleife und einer 2,4-µL Probennadel beträgt die Obergrenze für die Probenmenge, die noch vollständig injiziert werden kann, 6,4 µL. Wird ein höheres Injektionsvolumen gewählt, wird in Chromeleon eine Warnung ausgegeben. Das gewählte Injektionsvolumen wird zwar aufgezogen, aber die Probe wird nicht vollständig injiziert und die Genauigkeit ist etwas geringer.

Ein Beispiel für benutzerdefinierte Programme mit Mikroliter Pick-up finden Sie in der Hilfe zu Chromeleon.

Nachfolgend ist die Schaltsequenz beim Mikroliter-Pickup beschrieben:

Zu Beginn der Sequenz befindet sich das Ventil in der Position INJECT ( $\rightarrow$  Abb. 80). Der Vorstecher hat das Septum des Transportfläschchens durchstoßen und die Probennadel befindet sich im Fläschchen.

**I** Hinweis: Es wird empfohlen, bei Mikroliter-Pickup-Injektionen die mobile Phase als Transportflüssigkeit zu verwenden. So vermeiden Sie, dass die Transportflüssigkeit im Chromatogramm als separater Peak erscheint.



Abb. 80: Position INJECT

Bei der ersten Injektion nach einem Waschzyklus oder nachdem die Pufferschleife entleert wurde, zieht die Spritze die Transportflüssigkeit aus dem Transportfläschchen auf. Indem die Probennadel mit Transportflüssigkeit gefüllt wird ( $\rightarrow$  Abb. 81), wird die restliche Waschflüssigkeit, die sich eventuell noch in der Kapillare befindet, daraus entfernt.



Abb. 81: Aufziehen der Transportflüssigkeit

Die Probennadel verlässt das Transportfläschchen und fährt in das Probengefäß. Das Injektionsventil wird in die Position LOAD geschaltet ( $\rightarrow$  Abb. 82).



Abb. 82: Position LOAD

Das programmierte Injektionsvolumen wird aus dem Probengefäß aufgezogen ( $\rightarrow$  Abb. 83).



Abb. 83: Aufziehen des Probenvolumens

Die Probennadel fährt in das Transportfläschchen zurück ( $\rightarrow$  Abb. 84). Die Probenmenge wird mit der Transportflüssigkeit (mobile Phase) aus dem Transportfläschchen in die Probenschleife transportiert.



Abb. 84: Transportieren der Probe in die Probenschleife

Das Injektionsventil wird in die Position INJECT geschaltet ( $\rightarrow$  Abb. 85). Die Probenschleife wird Teil des Flussweges der mobilen Phase; die Probe wird auf die Säule transportiert und die Analysenzeit beginnt.



Abb. 85: Position INJECT

Die folgende Abbildung zeigt die Verteilung der Flüssigkeiten beim Mikroliter-Pickup:





**i** Hinweis: Wenn nach der Injektion ein Waschzyklus durchgeführt wird oder wenn der Inhalt der Pufferschleife in den Wasch-Port entleert wurde, wird die komplette Sequenz wiederholt, also auch Transportflüssigkeit zu Beginn der Sequenz aufgezogen. In allen anderen Fällen entfällt dieser Schritt.

# 8.4.4 Low Dispersion-Modus

Zur Optimierung des Injektionsprofils unterstützt der Autosampler für Full-Loop- und Partial-Loop-Injektionen den Low Dispersion-Modus. Die entsprechenden Kommandos und Eigenschaften dafür stehen in Chromeleon zur Verfügung.

Hinweis: Der Low Dispersion-Modus kann auch in benutzerdefinierten Programmen (UDP) verwendet werden.

Das Injektionsventil wird nach einer berechneten Zeit wieder in die Position LOAD zurückgeschaltet; die Rückflanke des Probenpfropfens wird abgeschnitten. Auf diese Weise können Sie das Injektionsprofil optimieren und damit auch die Reproduzierbarkeit verbessern.

Folgende Parameter sind wichtig:

#### • LowDispersionFlow:

Dieser Wert ist die Flussrate, mit der die Probe aus der Probenschleife heraus gespült wird.

**Hinweis:** Wird die Flussrate auf 0.00 µL/min gesetzt, wird eine Standardinjektion durchgeführt.

#### • LowDispersionFactor (L.D.-Faktor):

Über **LowDispersionFactor** legen Sie das Probenvolumen fest, das abgeschnitten wird, wenn das Injektionsventil zurückgeschaltet wird. Der Wert 1,0 entspricht dabei dem Volumen des Probenpfropfens. Ist der Wert höher (niedriger) als 1, wird das Injektionsventil zurückgeschaltet, wenn das Volumen größer (kleiner) ist als das Probenvolumen, das über die Probenschleife transportiert wurde. Der gültige Bereich beträgt 0,01 bis 100. Wird beispielsweise ein Wert von 1,2 gesetzt, wird das Injektionsventil zurückgeschaltet, wenn 120% des Probenpfropfens transportiert worden sind. Abb. 87 zeigt das Injektionsprofil bei unterschiedlichen Werten:



Abb. 87: Low Dispersion-Injektionsprofil

Mit folgender Formel wird die Zeit berechnet, nach der das Injektionsprofil in die Position LOAD zurückgeschaltet wird:

St = (f) (V1) (60)/F

wobei:  $S_t = Schaltzeit (Sek)$ 

f = L.D.-Faktor (Low Dispersion-Faktor)  $V_1 = Injektionsvolumen(\mu L)$  $F = LowDispersionFlow (\mu L/min)$ 

Ist beispielsweise das Injektionsvolumen  $V_1 = 1 \ \mu L$  und die Flussrate  $F = 5 \ \mu L$  und ist der L.D.-Faktor f = 1,1, beträgt die Schaltzeit  $S_t = 13,2$  Sek.

**I** Hinweise: Unter diesen Bedingungen wird die Probenschleife auch mit der Waschflüssigkeit gespült. Dies kann zu Injektionsproblemen führen, beispielsweise zu einer nicht einwandfreien Auflösung der Peaks. Wir empfehlen daher, als Waschflüssigkeit die mobile Phase (typischerweise die mobile Phase A) zu verwenden.

Die Schaltzeit ( $S_t$ ) muss mindestens 10 Sek. betragen. Liegt die berechnete Schaltzeit unter diesem Wert, wird sie auf 10 Sek. gesetzt.

# 9 Technische Daten

	WPS-3000(T)PL (RS)/ WPS-3000TBPL Nano/Cap	WPS-3000TBPL Analytisch/ WPS-3000T(B)FC	
Probenkapazität	Abhängig von der Drehtellerkonfiguration:		
Probenfläschchen	216 x 0,3 mL-Probenfläschchen, 120 x 1,1 mL-Probenfläschchen (konisch), 216 x 1,2 mL-Probenfläschchen, 120 x 1,8 mL- oder 2,0 mL- Probenfläschchen, 66 x 4 mL-Probenfläschchen, 30 x 10 mL- Probenfläschchen und/oder 120 x 0,5 mL- oder 1,5 mL-Eppendorf-Gefäße (offen, ohne Kappe) + 15 x 10 mL-Probenfläschchen		
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		
Injektionsmethoden	Full-Loop- und Partial-Loop-Injektionen (beide auch im Low-Dispersion-Modus) Mikroliter Pick-up Frei programmierbare, anwenderdefinierte Injektionsmethoden		
Injektionstechnik	Nadel in Nadel, programmierbares Waschen der Nadel		
Waschen der Nadel	Aktives Waschen der Probennadel von außen		
Verschleppung	< 0,02% für Koffein mit externem Waschen		
Erforderliche Probenmenge (Minimum)	1μL von 1μL (Mikroliter Pick-up)		
Injektionsvolumen Einstellbar: Empfohlen:	0,001 - 20 μL 20 nL - 20 μL	0,01 - 250 μL 0,1 - 250 μL	
Fraktioniermodi		nur WPS-3000T(B)FC: zeitabhängig, peakabhängig, manuelle und externe Trigger	
Präzision des Injektionsvolumens	RSD < 0,4% bei 1 μL-Full-Loop-Injektionen RSD < 1% bei 200 nL-Partial-Loop- Injektionen	RSD < 0,25% bei 5 μL-Full-Loop-Injektionen RSD < 0,3% bei 5μL/20 μL-Partial- Loop-Injektionen	
Linearität	Korr. koeff. > 0,9995, bei 100 nL- bis 500 nL- Partial-Loop-Injektionen, Koffein in Wasser	Korr. koeff. > 0,9999, RSD < 0,5 % bei 5 - 30 $\mu$ L (Koffein in Wasser), Offset < 2,0 $\mu$ L (500 $\mu$ L-Puffer- schleife), < 11,0 $\mu$ L (1000 $\mu$ L- Pufferschleife)	
Injektionsdauer	< 30 Sekunden bei 1 µL-Full-Loop-Injektionen	< 60 Sekunden bei 5 μL-Full-Loop-Injektionen < 90 Sekunden bei 5 μL-Partial-Loop-Injektionen	

	WPS-3000(T)PL (RS)/ WPS-3000TBPL Nano/Cap	WPS-3000TBPL Analytisch/ WPS-3000T(B)FC	
WPS-3000TPL (RS), TBPL und TBFC			
Probentemperierung Genauigkeit Probentemperatur	4 bis 45°C, bis zu 22°C unterhalb der $\pm 2$ °C bei einem Sollwert von 10 °C be °C und $\leq 50$ % relativer Luftfeuchtigke	Umgebungstemperatur ei Umgebungstemperaturen von≤ 25 eit	
Steuerung	Alle Funktionen softwaregesteuert übe Integrierter USB 1.1 Hub mit 3 freien	er USB 1.1 USB-Ports	
E/A-Schnittstellen	4 digitale Eingänge, 4 programmierbar	re Relaisausgänge	
Eingabe/Anzeige	LCD zur Anzeige von Systemparametern Standby-Schalter 3 LEDs (Power, Connected und Status) zur Statusangabe 4 Softkeys zum Betrieb während der Erstinstallation und zur Wartung		
Sicherheitsmerkmale	Leaksensor, automatische Mikrotiterplatten-/Fläschchenerkennung, Überlastungsschutz der Probennadel (Vorstecher), interne Überwachung aller mechanischen Operationen Autosampler mit Probenthermostatisierung: Überwachung der Funktion und Kühlleistung über 3 Temperatursensoren		
GLP	Unter Chromeleon: Automatische Gerätequalifikation(AutoQ <sup>TM</sup> -Routinen) und Überwachung der Wellness. Alle Systemparameter werden im Audit Trail von Chromeleon protokolliert.		
Medienberührte Teile			
Flusspfad Probe	PEEK, PAEK, PEEKsil, Fused Silica, Edelstahl	PEEK, PAEK, PCTFE Edelstahl (WPS-3000TFC)	
Flusspfad Waschflüssigkeit	PCTFE, ETFE, ECTFE, PTFE, Glas	PCTFE, ETFE, ECTFE, PTFE, Glas	
Leistungsaufnahme	100–120 V, 60 Hz; 200–240 V, 50 Hz; Automatische Spannungswahl max. 150 W (WPS-3000PL) bzw. max. 320 W (WPS-3000TPL/TFC)		
Emissionsschalldruckpegel	Typisch < 65 dB(A) in 1 m Abstand		
Umgebungsbedingungen	Verwendungsbereich:InnenraumTemperaturbereich:10 °C bis 35 °CLuftfeuchtigkeit:80% relative Feuchte, nicht kondensierendÜberspannungskat.:IIVerschmutzungsgrad:2		
Abmessungen	$36 \times 42 \times 51 \text{ cm} (\text{H} \times \text{B} \times \text{T})$		
Gewicht	WPS-3000PL (RS): 19 kg WPS-3000TPL (RS): 24 kg WPS-3000TBPL Nano/Cap: 24 kg	WPS-3000TBPL Analytisch: 24 kg WPS-3000T(B)FC: 25 kg	

Technische Daten: September 2013. Änderungen vorbehalten!

# 10 Zubehör, Ersatzteile und Verbrauchsmaterialien

Ersatzteile und Zubehör 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 Zubehör 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 180).

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-Vertriebsorganisation für Dionex HPLC-Produkte.

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000PL RS, WPS-3000TPL RS		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	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 (250V 6,3 x 32 mm)	Enthalten in 6820.0026	1
Innensechskantschlüssel, Größe 9/64"	6000.0053	1
Gabelschlüssel, Größe 1/4 x 5/16"	6000.0051	1
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011	1
Probennadel (2,4 µL, Fused Silica, nanoViper; 100 µm ID)	6820.3010	1
Probenschleife (20 µL, nanoViper)	6826.2420	1
Probenhalter für 40 Probenfläschchen, A.D. < 12 mm und Standard-Mikrotiterplatten (Höhe 12 - 24 mm)	6820.4070	3
Transportfläschchen (10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Waschflasche WPS	Enthalten in 6820.4075	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102	1

Bezeichnung	BestNr.	Menge im Zubehör
Polypropylen-Fläschchen (250 µL)	Enthalten in 6820.0029	10
Deckel für 250 µL Polypropylen-Fläschchen	Enthalten in 6820.0028	10
USB-Kabel Typ A auf Typ B (USB 2.0, 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000PL und WPS-3000TPL Nano/Cap mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	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 (250V 6,3 x 32 mm)	Enthalten in 6820.0026	1
Innensechskantschlüssel, Größe 9/64"	6000.0053	1
Gabelschlüssel, Größe 1/4 x 5/16"	6000.0051	1
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011	1
Probennadel (2,4 µL, Fused Silica, nanoViper; 100 µm ID)	6820.3010	1
Probenschleife (20 µL, nanoViper)	6826.2420	1
Probenhalter für 40 Probenfläschchen, A.D. < 12 mm und Standard-Mikrotiterplatten (Höhe 12 - 24 mm)	6820.4070	3
Transportfläschchen (10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Waschflasche WPS	Enthalten in 6820.4075	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102	1
Polypropylen-Fläschchen (250 µL)	Enthalten in 6820.0029	10
Deckel für 250 µL Polypropylen-Fläschchen	Enthalten in 6820.0028	10
USB-Kabel Typ A auf Typ B (USB 2.0, 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000TBPL Nano/Cap (biokompatibler Autosampler) mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	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 (250V 6,3 x 32 mm)	Enthalten in 6820.0026	1
Innensechskantschlüssel, Größe 9/64"	6000.0053	1
Gabelschlüssel, Größe 1/4 x 5/16"	6000.0051	1
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011	1
Probennadel (2,4 µL, Fused Silica, nanoViper; 100 µm ID)	6820.3010	1
Probenschleife (20 µL, PEEKsil)	6821.0018	1
Probenhalter (für 40 Probenfläschchen, A.D. < 12 mm)	6820.4070	3
Transportfläschchen (10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Waschflasche WPS	Enthalten in 6820.4075	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102	1
Polypropylen-Fläschchen (250 µL)	Enthalten in 6820.0029	10
Deckel für 250 µL Polypropylen-Fläschchen	Enthalten in 6820.0028	10
USB-Kabel Typ A auf Typ B (USB 2.0, 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000TBPL Analytisch (biokompatibler Autosampler) mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	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 (250V 6,3 x 32 mm)	Enthalten in 6820.0026	1
Werkzeug zur Fittingvormontage	6000.0065	1
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011	1
Kapillare (PEEK, 1/16" AD x 0,25 mm ID) WPS - TCC	6251.6001	2 m
Einteiliges Fitting (PEEK, 1/16", 15 mm)	6266.0024	2
Probennadel (15 µL, PEEK, Viper; 250 µm ID)	6820.3025	1
Probenschleife (250 µL, PEEK)	6823.0020	1
Finger-tight-Fittingschraube (PEEK, 1/16", mit Ferrule) zum Anschluss von Kapillaren am Injektionsventil	6823.0010	2
Pufferschleife (1000 µL, PEEK/PTFE)	6821.0022	1
Spritze (250 µl)	6822.0003	1
Probenhalter (für 40 Probenfläschchen, A.D. < 12 mm)	6820.4070	3
Transportfläschchen (10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Probenfläschchen (1,8 mL, braun)	Enthalten in 6000.0072	10
Probenfläschchendeckel mit Septum für 1,8 mL-Fläschchen	Enthalten in 6000.0071	10
Probenfläschchen (2 mL, braun)	Enthalten in 6000.0060	10
Probenfläschchendeckel (weiß) für 2 mL-Probenfläschchen	Enthalten in 6000.0057	10
Probenfläschchenseptum (Silikon, rot) für 2 mL-Fläschchen	Enthalten in 6000.0058	10
Waschflasche WPS	Enthalten in 6820.4075	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102	1
USB-Kabel Typ A auf Typ B (USB 2.0, 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000TFC mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	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 (250V 6,3 x 32 mm)	Enthalten in 6820.0026	1
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011	1
Probennadel (15µL, Fused Silica, nanoViper; 250 µm ID)	6820.3115	1
Probenschleife (250 µL, Stahl)	6824.0020	1
Pufferschleife (1000 µL, Stahl/PTFE)	6820.0056	1
Probenhalter für 40 Probenfläschchen, A.D. < 12 mm und Standard-Mikrotiterplatten (Höhe 12 - 24 mm)	6820.4070	3
Transportfläschchen (10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Waschflasche WPS	Enthalten in 6820.4075	1
Probenfläschchen (1,8 mL, braun)	Enthalten in 6000.0072	10
Probenfläschchendeckel mit Septum für 1,8 mL-Fläschchen	Enthalten in 6000.0071	10
Probenfläschchen (2 mL, braun)	Enthalten in 6000.0060	10
Probenfläschchendeckel (weiß) für 2 mL-Probenfläschchen	Enthalten in 6000.0057	10
Probenfläschchenseptum (Silikon, rot) für 2 mL-Fläschchen	Enthalten in 6000.0058	10
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Deep-Well-Plate (34 - 46 mm Höhe, 96 Kavitäten)	6820.4101	1
Probenhalter für Deep-Well-Plates (34 - 46 mm Höhe)	6820.4079	1
Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102	2
USB-Kabel Typ A auf Typ B (USB 2.0, 1 m)	6035.9035	1

Bezeichnung	BestNr.	Menge im Zubehör
Zubehör für WPS-3000TBFC mit:		
Online Degas Wash-Kit (Enthält alle Schläuche und Fittinge, die für den Anschluss der Waschflüssigkeit zur Online-Entgasung benötigt werden.)	6820.2450	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 (250V 6,3 x 32 mm)	Enthalten in 6820.0026	1
Werkzeug zur Fittingvormontage	6000.0065	1
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011	1
Kapillare (PEEK, 1/16" AD x 0,25 mm ID) für WPS - TCC bzw. Pumpe - WPS	6251.6001	2 m
Einteiliges Fitting (PEEK, 1/16", 15 mm)	6266.0024	2
Probennadel (15 µL, PEEK, Viper, 250 µm ID)	6820.3025	1
Probenschleife (250 µL, PEEK)	6823.0020	1
Fingertight-Fitting (PEEK, 1/16", mit Ferrule) zum Anschluss von Kapillaren am Injektionsventil	6823.0010	2
Pufferschleife (1000 µL, PEEK/PTFE)	6821.0022	1
Probenhalter (für 40 Probenfläschchen, A.D. < 12 mm)	6820.4070	3
Transportfläschchen (10 mL) mit Kappe und Septum	Enthalten in 6820.0023	16
Probenfläschchen (1,8 mL, braun)	Enthalten in 6000.0072	10
Probenfläschchendeckel mit Septum für 1,8 mL-Fläschchen	Enthalten in 6000.0071	10
Waschflasche WPS	Enthalten in 6820.4075	1
Standard-Mikrotiterplatte (96 Kavitäten)	6820.4100	1
Deep-Well-Plate (34 - 46 mm Höhe, 96 Kavitäten)	6820.4101	1
Probenhalter für Deep-Well-Plates (34 - 46 mm Höhe)	6820.4079	1
Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102	2
USB-Kabel Typ A auf Typ B (USB 2.0, 1 m)	6035.9035	1

# 10.2 Optionales Zubehör

# 10.2.1 WPS-3000(T)PL (RS)

Optionales Zubehör	BestNr.	Bemerkungen
Automated Off-Line 2DLC Solution	6820.0050	Micro Fraction Collection Option plus Chromeleon Timebase Class 1 und Extended Fraction Collection License
Drainage-Kit für UltiMate 3000 Systeme	6040.0005	Das Kit enthält alle erforderlichen Komponenten für den Anschluss der Systemdrainage sowie eine detaillierte Installationsanleitung.
Extended Fraction Collection License	5960.0039	Chromeleon-Zusatzlizenz für die erweiterte Fraktionssammlung (→ Seite 141) Diese Lizenz schließt die Fraction Collection License ein.
Fraction Collection License	5960.0038	Chromeleon-Zusatzlizenz mit Grundfunktionen für das Sammeln von Fraktionen (→ Seite 141)
Frontabdeckung (transparent)	6820.1427	Verschließt den Probenraum zum Schutz gegen Staub oder andere Partikel.
Micro Fraction Collection Option	6820.0051	Erweiterung (8-Port-Ventil) zum Sammeln von Fraktionen (ohne Software-Lizenz).
PAEK Modification Kit	Wenden Sie sich an den Service.	Zur Umrüstung der nicht- biokompatiblen Fluidik eines WPS-3000(T)PL RS Autosamplers zu einem biokompatiblen Autosampler
Optionales Zubehör für biokompatible Anwendungen mit WPS-3000(T)PL (RS) Nano/Kap:		
<ul> <li>Probenschleifen (PEEK)</li> <li>5 μL</li> <li>10 μL</li> <li>20 μL</li> <li>50 μL, für analytische Anwendungen</li> <li>125 μL, für analytische Anwendungen</li> </ul>	6823.0016 6823.0017 6823.0018 6823.0019 6821.0032	
Spritzen-Bio-Kit, enthält Spritze, 250 μL Nadel, 15 μL, Fused Silica, nanoViper Pufferschleife, 500 μL, PEEK/PTFE Probenschleife, 125 μL, PEEK	6821.0031	
<i>Pufferschleifen, PEEK/PTFE</i> 50 μL 500 μL	6821.0019 6821.0020	

Optionales Zubehör	BestNr.	Bemerkungen
Probennadeln		
3,5 $\mu$ L, PEEK, Viper; 127 $\mu$ m ID 15 $\mu$ L, PEEK, Viper; 250 $\mu$ m ID; für analytische Zwecke	6820.3023 6820.3025	
für analytische Zwecke	6820.3115	
Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 1 μL 5 μL 10 μL 20 μL 50 μL 125 μL, für analytische Anwendungen	6826.2401 6826.2405 6826.2410 6826.2420 6826.2450 6826.2412	Vergewissern Sie sich, dass auch eine geeignete Pufferschleife und Spritze installiert ist (→ Seite 16).
<i>Pufferschleife, SST/PTFE</i> 500 μL	6820.0020	Vergewissern Sie sich, dass auch eine geeignete Spritze und Probenschleife installiert ist (→ Seite 16).
Signalkabel (6-pol. Mini-DIN)	6000.1004	Anschluss von externen Geräten.
Spritzen 100 μL 250 μL, für analytische Anwendungen	6822.0002 6822.0003	Vergewissern Sie sich, dass auch die geeigneten Puffer- und Probenschleifen installiert sind ( $\rightarrow$ Seite 16).
Spritzenkit (250 µL), enthält Spritze, 250 µL Nadel, 15 µL, Fused Silica, nanoViper, Pufferschleife, 500 µL, SST/PTFE Probenschleife, 125 µL, nanoViper	6820.0031	Komplett-Kit zum Aufrüsten auf eine 250 µL-Spritze
## 10.2.2 WPS-3000TBPL Nano/Cap

Optionales Zubehör	BestNr.	Bemerkungen
Automated Off-Line 2DLC Solution (biokompatibel)	6821.0050	Micro Fraction Collection Option plus Chromeleon Timebase Class 1 und Extended Fraction Collection License
Drainage-Kit für UltiMate 3000 Systeme	6040.0005	Das Kit enthält alle erforderlichen Komponenten für den Anschluss der Systemdrainage sowie eine detaillierte Installationsanleitung.
Extended Fraction Collection License	5960.0039	Chromeleon-Zusatzlizenz für die erweiterte Fraktionssammlung (→ Seite 141) Diese Lizenz schließt die Fraction Collection License ein.
Fraction Collection License	5960.0038	Chromeleon-Zusatzlizenz mit Grundfunktionen für das Sammeln von Fraktionen (→ Seite 141)
Frontabdeckung (transparent)	6820.1427	Verschließt den Probenraum zum Schutz gegen Staub oder andere Partikel.
Micro Fraction Collection Option (biokompatibel)	6821.0051	Erweiterung (8-Port-Ventil) zum Sammeln von Fraktionen (ohne Software-Lizenz).
<i>Probennadeln</i> 15 μL, PEEK, Viper; 250 μm ID; für analytische Zwecke 15 μL, Fused Silica, nanoViper; 250 μm ID; für analytische Zwecke	6820.3025 6820.3115	
Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 1 μL 5 μL 10 μL 20 μL 50 μL, für analytische Anwendungen 125 μL, für analytische Anwendungen Probenschleifen, PEEK 5 μL 10 μL 20 μL 50 μL 125 μL	6826.2401 6826.2405 6826.2410 6826.2420 6826.2450 6826.2412 6823.0016 6823.0017 6823.0018 6823.0019 6821.0032	Vergewissern Sie sich, dass auch eine geeignete Pufferschleife und Spritze installiert ist (→ Seite 16).
<i>Pufferschleife, PEEK/PTFE</i> 500 μL	6821.0020	Vergewissern Sie sich, dass auch die geeigneten Spritzen und Probenschleifen installiert sind (→ Seite 16).
Signalkabel (6-pol. Mini-DIN)	6000.1004	Anschluss von externen Geräten.

Optionales Zubehör	BestNr.	Bemerkungen
Spritzen 100 μL 250 μL, für analytische Anwendungen	6822.0002 6822.0003	Vergewissern Sie sich, dass auch die geeigneten Puffer- und Probenschleifen installiert sind ( $\rightarrow$ Seite 16).
Spritzen-Bio-Kit, enthält Spritze, 250 µL Nadel, 15 µL, Fused Silica, nanoViper Pufferschleife, 500 µL, PEEK/PTFE Probenschleife, 125 µL, PEEK	6821.0031	Komplett-Kit (biokompatibel) zum Aufrüsten auf eine 250 µL-Spritze

10.2.3 WP	S-3000TBPL Analytisch
-----------	-----------------------

Optionales Zubehör	BestNr.	Bemerkungen
Drainage-Kit für UltiMate 3000 Systeme	6040.0005	Das Kit enthält alle erforderlichen Komponenten für den Anschluss der Systemdrainage sowie eine detaillierte Installationsanleitung.
Frontabdeckung (transparent)	6820.1427	Verschließt den Probenraum zum Schutz gegen Staub oder andere Partikel.
Probennadeln		
3,5 μL, PEEK, Viper; 127 μm ID 15 μL, Fused Silica, nanoViper; 250 μm ID	6820.3023	
	6820.3115	
Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 50 μL 125 μL Probenschleifen, PEEK 5 μL 10 μL 20 μL 50 μL 250 μL	6826.2450 6826.2412 6823.0016 6823.0017 6823.0018 6823.0019 6821.0032 6823.0020	Vergewissern Sie sich, dass auch eine geeignete Pufferschleife und Spritze installiert sind (→ Seite 16). Die Probenschleifen mit 5, 10 und 20 µL sind für Full-Loop-Injektionen mit diesen Volumina vorgesehen.
<i>Pufferschleife, PEEK/PTFE</i> 1000 μL	6821.0022	Vergewissern Sie sich, dass auch eine geeignete Spritze und Probenschleife installiert ist ( $\rightarrow$ Seite 16).
Signalkabel (6-pol. Mini-DIN)	6000.1004	Anschluss von externen Geräten.
Spritze, 250 µL	6822.0003	Vergewissern Sie sich, dass auch die geeigneten Puffer- und Probenschleifen installiert sind ( $\rightarrow$ Seite 16).

## 10.2.4 WPS-3000TFC

Optionales Zubehör	BestNr.	Bemerkungen
Drainage-Kit für UltiMate 3000 Systeme	6040.0005	Das Kit enthält alle erforderlichen Komponenten für den Anschluss der Systemdrainage sowie eine detaillierte Installationsanleitung.
Extended Fraction Collection License	5960.0039	Chromeleon-Zusatzlizenz für die erweiterte Fraktionssammlung (→ Seite 141) Diese Lizenz schließt die Fraction Collection License ein.
Fraction Collection License	5960.0038	Chromeleon-Zusatzlizenz mit Grundfunktionen für das Sammeln von Fraktionen (→ Seite 141)
Frontabdeckung (transparent)	6820.1427	Verschließt den Probenraum zum Schutz gegen Staub oder andere Partikel.
<ul> <li>Kit zur Optimierung für Nano/Kap- Anwendungen mit</li> <li>2.4 μL-Probennadel (Fused Silica, nanoViper)</li> <li>25 μL-Spritze</li> <li>50 μL-Pufferschleife</li> <li>1 μL-Probenschleife, nanoViper</li> <li>1 μL-Bridge (PEEKsil)</li> </ul> Optionales Zubehör für Nano/Kap- Anwendungen: Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 1μL 5 μL 10 μL 20 μL Spritze 25 μL Pufferschleife, SST 50 μL Probennadeln 2,4 μL, Fused Silica, nanoViper; 100 μm ID	6824.0030 6826.2401 6826.2405 6826.2410 6826.2420 6822.0001 6820.0019 6820.3010	Installieren Sie die Komponenten aus dem Umrüstkit wie in den jeweiligen Servicekapiteln beschrieben (→ Kapitel 7, ab Seite 111 und 8.2.1.6, Seite 144)
3,5 μL, PEEK, Viper; 127 μm ID         Normalphasen-Kit, mit         - 15 μL-Probennadel, SST, Viper         - 6,4 μL-Bridge, SST, Viper         - Innensechskantschlüssel (2 mm)         - 1 Unterlegscheibe         - 1 Sicherungsscheibe	6820.0060	Für Normalphasen-Anwendungen mit einem WPS-3000TFC Autosampler. Das Kit enthält alle erforderlichen Komponente, um die Nadel, das Injektionsventil und das Fraktionierventil für Normalphasen- Anwendungen zu erden.

Optionales Zubehör	BestNr.	Bemerkungen
Probennadel		
15 μL, PEEK, Viper; 250 μm ID	6820.3025	
Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 50 μL 125 μL Probenschleife, SST 250 μL	6826.2450 6826.2412 6824.0020	Vergewissern Sie sich, dass auch eine geeignete Pufferschleife und Spritze installiert sind (→ Seite 16).
<i>Pufferschleife, SST/PTFE</i> 1000 μL	6820.0056	Vergewissern Sie sich, dass auch eine geeignete Probenschleife und Spritze installiert sind ( $\rightarrow$ Seite 16).
Signalkabel (6-pol. Mini-DIN)	6000.1004	Anschluss von externen Geräten.
Spritze, 100 µL	6822.0002	Vergewissern Sie sich, dass auch die geeigneten Puffer- und Probenschleifen installiert sind ( $\rightarrow$ Seite 16).

## 10.2.5 WPS-3000TBFC

Optionales Zubehör	BestNr.	Bemerkungen
Drainage-Kit für UltiMate 3000 Systeme	6040.0005	Das Kit enthält alle erforderlichen Komponenten für den Anschluss der Systemdrainage sowie eine detaillierte Installationsanleitung.
Extended Fraction Collection License	5960.0039	Chromeleon-Zusatzlizenz für die erweiterte Fraktionssammlung (→ Seite 141) Diese Lizenz schließt die Fraction Collection License ein.
Fraction Collection License	5960.0038	Chromeleon-Zusatzlizenz mit Grundfunktionen für das Sammeln von Fraktionen (→ Seite 141)
Frontabdeckung (transparent)	6820.1427	Verschließt den Probenraum zum Schutz gegen Staub oder andere Partikel.
<ul> <li>Nano/Kap-Umrüstkit, mit</li> <li>2,4 μL-Probennadel (Fused Silica, nanoViper)</li> <li>25 μL-Spritze</li> <li>50 μL-Pufferschleife, PEEK/PTFE</li> <li>1 μL-Probenschleife, nanoViper (Fused Silica/PEEK sheathed)</li> <li>1 μL-Bridge (PEEKsil, biokomp.)</li> </ul>	6825.0030	Installieren Sie die Komponenten aus dem Umrüstkit wie in den jeweiligen Servicekapiteln beschrieben (→ Kapitel 7, ab Seite 111 und 8.2.1.6, Seite 144)
Anwendungen: Probenschleifen, nanoViper (Fused Silica/PEEK sheathed)		
1 μL 5 μL 10 μL 20 μL	6826.2401 6826.2405 6826.2410 6826.2420	
Spritze 25 μL	6822.0001	
Pufferschleife, PEEK/PIFE 50 μL	6821.0019	
2,4 μL, Fused Silica, nanoViper; 100 μm ID 3,5 μL, PEEK, Viper; 127 μm ID	6820.3010 6820.3023	
Probennadeln		
3,5 μL, PEEK, Viper; 127 μm ID 15 μL, PEEK, Viper; 250 μm ID	6820.3023 6820.3025	

Optionales Zubehör	BestNr.	Bemerkungen
Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 50 μL 125 μL Probenschleifen, PEEK 5 μL 10 μL 20 μL 50 μL 125 μL 250 μL	6826.2450 6826.2412 6823.0016 6823.0017 6823.0018 6823.0019 6821.0032 6823.0020	Vergewissern Sie sich, dass auch eine geeignete Pufferschleife und Spritze installiert sind (→ Seite 16).
<i>Pufferschleife, PEEK/PTFE</i> 1000 μL	6821.0022	Vergewissern Sie sich, dass auch eine geeignete Probenschleife und Spritze installiert sind ( $\rightarrow$ Seite 16).
Signalkabel (6-pol. Mini-DIN)	6000.1004	Anschluss von externen Geräten.
Spritze, 100 µL	6822.0002	Vergewissern Sie sich, dass auch die geeigneten Puffer- und Probenschleifen installiert sind ( $\rightarrow$ Seite 16).

# **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-Vertriebsorganisation für Dionex HPLC-Produkte.

Bezeichnung	BestNr.
Adapter für 2 mL-Probenfläschchen in 10 mL-Transportfläschchenhalter	6820.4092
Adapter für Low-Well-PCR-Plates mit einer Höhe von 8 - 12 mm	6820.4088
Adapter für Niederhalter bei Verwendung von 24-Deep-Well-Plates	6820.2402
Bridge (PEEK, Volumen: 6,2 µL) für WPS-3000TFC WPS-3000TBFC	6824.0016 6825.0016
Bridge (PEEKsil, Volumen: 1 µL) für WPS-3000TFC, optimiert für Nano/Kap-Anwendungen WPS-3000TBFC, optimiert für Nano/Kap-Anwendungen	6824.0015 6825.0015
Bridge für WPS-3000TFC für Normalphasen-Anwendungen (Kapillare, Viper, SST, Volumen 6,4 $\mu$ L, 0,18 x 250 mm (I.D. x L)) Enthalten im Normalphasen-Kit (BestNr. 6820.0060). Informationen zum Kit finden Sie in Kapitel 10.2.4 ( $\rightarrow$ Seite 176).	6040.2385
Dichtung Spritze/Injektionsventil (5 Stück)	6822.0009
Drainage-Kit für UltiMate 3000 Systeme	6040.0005
Einteiliges Fitting (PEEK, 1/16", 15 mm)	6266.0024
Finger-tight-Fittingschraube (PEEK, 1/16") zum Anschluss der Kapillaren am Injektionsventil des WPS-3000TBPL Analytisch	6823.0010
Frontabdeckung (transparent)	6820.1427
Gabelschlüssel, Größe 1/4 x 5/16"	6000.0051
Innensechskantschlüssel, Größe 9/64"	6000.0053
Kapillare (kurz) zwischen Ladepumpe und Autosampler, biokompatible Geräte (130 µm ID x 60 cm, PEEK), einschl. geeigneter Fittingverbindungen	6721.0031
Kapillare (kurz) zwischen Ladepumpe und Autosampler, Standardgeräte (130 µm ID x 60 cm, PEEK), einschl. geeigneter Fittingverbindungen	6720.0031
Kapillare (lang) zwischen Ladepumpe und Autosampler, biokompatible Geräte (130 µm ID x 75 cm, PEEK), einschl. geeigneter Fittingverbindungen	6721.0032
Kapillare (lang) zwischen Ladepumpe und Autosampler, Standardgeräte (130 µm ID x 75 cm, PEEK), einschl. geeigneter Fittingverbindungen	6720.0032
Kapillare (PEEK, 1/16" AD x 0,25 mm ID) zwischen WPS-3000TBPL Analytisch/WPS-3000TBFC und Säulenthermostat bzw. zwischen Pumpe und WPS-3000TBPL Analytisch	6251.6001
Kapillare zwischen Ladepumpe und Autosampler (130 µm ID x 45 cm, PEEK) einschließlich geeigneter Fittingverbindungen, Standardgeräte	6720.0057

Bezeichnung	BestNr.
Kapillare zwischen Ladepumpe und Autosampler (130 µm ID x 45 cm, PEEK) einschließlich geeigneter Fittingverbindungen, biokompatible Geräte	6721.0057
Kapillare zwischen Pumpe und WPS-3000TFC (Viper, SST, 0,18 mm x 550 mm ID x L)	6040.2355
Kapillare zwischen WPS-3000TFC und TCC-3000 (Viper, SST, 0,13 mm x 350 mm ID x L)	6040.2335
Kapillar-Kupplung (1/16", totvolumenfrei)	6720.0039
Menüstift	6300.0100
Mikrotiterplatte (Deep-Well-Plate mit 384 Kavitäten)	Auf Anfrage
Mikrotiterplatte (Deep-Well-Plate mit 96 Kavitäten)	6820.4101
Mikrotiterplatte (Standard-Mikrotiterplatte mit 384 Kavitäten)	6820.4110
Mikrotiterplatte (Standard-Mikrotiterplatte mit 96 Kavitäten)	6820.4100
Mikrotiterplatte-Verschlussmatte für Mikrotiterplatte und Deep-Well-Plate mit 96 Kavitäten	6820.4102
Mikrotiterplatte-Verschlussmatte (für Standard-Mikrotiterplatten mit 384 Kavitäten)	6820.4112
Mutter (1/16", Standard, SS316, 10 Stück)	754.ZN1S6
Mutter (1/16", Standard, SS316, 10 Stück)	6720.0018
Netzkabel, Australien, China	6000.1060
Netzkabel, Dänemark	6000.1070
Netzkabel, EU	6000.1000
Netzkabel, Indien/SA	6000.1090
Netzkabel, Italien	6000.1040
Netzkabel, Japan	6000.1050
Netzkabel, Schweiz	6000.1030
Netzkabel, UK	6000.1020
Netzkabel, US	6000.1001
Normalphasen-Kit für WPS-3000TFC Autosampler Für Normalphasen-Anwendungen mit einem WPS-3000TFC Autosampler. Das Kit enthält alle erforderlichen Komponente, um die Nadel, das Injektionsventil und das Fraktionierventil für Normalphasen-Anwendungen zu erden. Informationen zum Inhalt des Kits finden Sie auf Seite 176.	6820.0060
Probenfläschchen (1,8 mL, braun, 100 Stück)	6000.0072
Probenfläschchen (4 mL; braun; 100 Stück)	6000.0074
Probenfläschchen, zylindrisch 1,2 mL (500 Stück)	6000.0062
Probenfläschchen, zylindrisch 2 mL (500 Stück), braun	6000.0060
Probenfläschchendeckel (Schraubdeckel) für 4-mL-Probenfläschchen (100 Stück)	6000.0073
Probenfläschchendeckel (weiß) für 2 mL-Probenfläschchen (500 Stück)	6000.0057

Bezeichnung	BestNr.
Probenfläschchendeckel für 1,2-mL-Fläschchen (Crimpkappe; 1000 Stück)	6000.0064
Probenfläschchendeckel mit Septum für 1,1 mL-Fläschchen (Bördelkappe mit Silikon- Septum, 1000 Stück)	6000.0076
Probenfläschchendeckel mit Septum für 1,8 mL-Fläschchen (Bördelkappe; 100 Stück)	6000.0071
Probenfläschchendeckel mit Septum für 1.2 mL-Fläschchen (Bördelkappe mit Silikon/PTFE-Septum, geschlitzt, 500 Stück)	6000.0061
Probenfläschchenseptum (Silikon, rot) für 2 mL-Fläschchen (500 Stück)	6000.0058
Probenfläschchenseptum (Silikon/PTFE, 500 Stück) für 4-mL-Probenfläschchen	6000.0075
Probenhalter für 10 zylindrische 10 mL Probenfläschchen	6820.4086
Probenhalter für 22 zylindrische 4 mL Probenfläschchen	6820.4084
Probenhalter für 40 Eppendorf-Gefäße 0,5 mL	6820.4096
Probenhalter für 40 Eppendorf-Gefäße 1,5 mL	6820.4094
Probenhalter für 40 konische 1.1 mL-Probenfläschchen	6820.4087
Probenhalter für 40 Probenfläschchen, A.D. < 12 mm	6820.4070
Probenhalter für 72 Mikrodialyse-Fläschchen	6820.4097
Probenhalter für 72 zylindrische 0,3 mL Probenfläschchen	6820.4091
Probenhalter für 72 zylindrische 1,2 mL Probenfläschchen	6820.4090
Probenhalter für Deep-Well-Plates 20 - 32 mm Höhe	6820.4089
Probenhalter für Deep-Well-Plates 30 - 36 mm Höhe	6820.4083
Probenhalter für Deep-Well-Plates 34 - 46 mm Höhe	6820.4079
<ul> <li>Probennadel für WPS-3000(T)PL (RS) Nano/Kap</li> <li>2,4 μL, Fused Silica, nanoViper; 100 μm ID</li> <li>3,5 μL, PEEK, Viper; 127 μm ID</li> <li>15 μL, PEEK, Viper; 250 μm ID; für analytische Zwecke</li> <li>15 μL, Fused Silica, nanoViper; 250 μm ID; für analytische Zwecke</li> </ul>	6820.3010 6820.3023 6820.3025 6820.3115
Probennadel für WPS-3000T(B)FC, optimiert für Nano/Kap-Anwendungen	(000 0010
2,4 μL, Fused Silica, nanoViper; 100 μm ID 3,5 μL, PEEK, Viper; 127 μm ID	6820.3010 6820.3023
Probennadel für WPS-3000TBFC 15 μL, PEEK, Viper; 250 μm ID 15 μL, Fused Silica, nanoViper; 250 μm ID	6820.3025 6820.3115
Probennadel für WPS-3000TBPL Analytisch 3,5 μL, PEEK, Viper; 127 μm ID 15 μL, PEEK, Viper; 250 μm ID 15 μL, Fused Silica, nanoViper; 250 μm ID	6820.3023 6820.3025 6820.3115

Bezeichnung	BestNr.
Probennadel für WPS-3000TBPL Nano/Kap 2,4 μL, Fused Silica, nanoViper; 100 μm ID 15 μL, PEEK, Viper; 250 μm ID; für analytische Zwecke 15 μL, Fused Silica, nanoViper; 250 μm ID; für analytische Zwecke	6820.3010 6820.3025 6820.3115
<ul> <li>Probennadel für WPS-3000TFC</li> <li>15 μL, PEEK, Viper; 250 μm ID</li> <li>15 μL, Fused Silica, nanoViper; 250 μm ID</li> <li>15 μL, SST, Viper; 200 μm ID, für Normalphasen-Anwendungen</li> <li>Enthalten im Normalphasen-Kit (BestNr. 6820.0060). Informationen zum Kit finden</li> <li>Sie in Kapitel 10.2.4 (→ Seite 176).</li> </ul>	6820.3025 6820.3115 
Probenschleife für WPS-3000(T)PL (RS) Nano/Cap, nanoViper (Fused Silica/PEEK sheathed) 1 μL 5 μL 10 μL 20 μL 50 μL 125 μL, für analytische Anwendungen	6826.2401 6826.2405 6826.2410 6826.2420 6826.2450 6826.2412
<ul> <li>Probenschleife für WPS-3000(T)PL (RS) Nano/Cap, umgerüstet mit PAEK Upgrade Kit</li> <li>5 μL, PEEK</li> <li>10 μL, PEEK</li> <li>20 μL, PEEK</li> <li>50 μL, PEEK, für analytische Anwendungen</li> <li>125 μL, PEEK, für analytische Anwendungen</li> </ul>	6823.0016 6823.0017 6823.0018 6823.0019 6821.0032
Probenschleife für WPS-3000TBFC Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 50 μL 125 μL Probenschleifen, PEEK 5 μL 10 μL 20 μL 50 μL 125 μL 250 μL	6826.2450 6826.2412 6823.0016 6823.0017 6823.0018 6823.0019 6821.0032 6823.0020
Probenschleife für WPS-3000TBPL Analytisch Probenschleifen, nanoViper (Fused Silica/PEEK sheathed) 50 μL 125 μL Probenschleifen, PEEK 5 μL 10 μL 20 μL 50 μL 125 μL 250 μL	6826.2450 6826.2412 6823.0016 6823.0017 6823.0018 6823.0019 6821.0032 6823.0020

Bezeichnung	BestNr.
Probenschleife für WPS-3000TBPL Nano/Cap Probenschleifen, nanoViper (Eused Silieg/PEEK sheathed)	
<ul> <li>Probenschleijen, hanoviper (Fused Stitca/PEEK sheathed)</li> <li>1 μL</li> <li>5 μL</li> <li>10 μL</li> <li>20 μL</li> <li>50 μL, für analytische Anwendungen</li> <li>125 μL, für analytische Anwendungen</li> </ul>	6826.2401 6826.2405 6826.2410 6826.2420 6826.2450 6826.2412
Probenschleifen, PEEK         5 μL         10 μL         20 μL         50 μL         125 μL	6823.0016 6823.0017 6823.0018 6823.0019 6821.0032
Probenschleife für WPS-3000TFC	
50 μL 125 μL	6826.2450 6826.2412
Probenschleifen, SST 50 μL 250 μL	6824.0019 6824.0020
Probenschleife für WPS-3000TFC und WPS-3000TBFC, optimiert für Nano/Kap- Anwendungen, nanoViper (Fused Silica/PEEK sheathed)	
1 μL 5 μL 10 μL 20 μL	6826.2401 6826.2405 6826.2410 6826.2420
Pufferschleife, PEEK/PTFE, für WPS-3000(T)PL (RS) Nano/Cap, umgerüstet mit PAEK Umrüst-Kit	
50 μL 500 μL	6821.0019 6821.0020
Pufferschleife, PEEK/PTFE, für WPS-3000TBFC 500 μL 1000 μL	6821.0020 6821.0022
Pufferschleife, PEEK/PTFE, für WPS-3000TBFC, optimiert für Nano/Kap- Anwendungen	
50 µL	6821.0019
Pufferschleife, PEEK/PTFE, für WPS-3000TBPL Analytisch 500 μL 1000 μL	6821.0020 6821.0022
Pufferschleife, PEEK/PTFE, für WPS-3000TBPL Nano/Cap	
50 μL 500 μL, für analytische Anwendungen	6821.0019 6821.0020

Bezeichnung	BestNr.
Pufferschleife, SST/PTFE, für WPS-3000(T)PL (RS) Nano/Cap 50 μL 500 μL	6820.0019 6820.0020
Pufferschleife, SST/PTFE, für WPS-3000TFC, optimiert für Nano/Kap-Anwendungen 50 μL	6820.0019
Pufferschleifen, SST/PTFE, für WPS-3000TFC 500 μL 1000 μL	6820.0020 6820.0056
Ringschlüssel für Vorstecher- und Fittingmontage	6000.0052
Rotor für Ventile mit 2-Positionen und 6 Ports Standardventil 900 bar-Ventil biokompatibles Ventil	6820.0014 6826.0013 6820.0035
Rotor für Ventile mit 2-Positionen und 8 Ports (Micro Fraction Collection Option) Standardventil (Edelstahl) biokompatibles Ventil	6820.0054 6821.0054
Schlauch zwischen Waschflasche und Spritzenventil/Abfall (ECTFE, 0,75 mm ID, 600 mm lang)	6820.0011
Schneidring (1/16", SST316, 10 Stück)	6720.0017
Schneidring- und Fittingkit (SR/FS-7), bestehend aus: 6 Fittingschrauben FS-7, 1/16", Edelstahl 6 Schneidringen SR-7, 1/16", Edelstahl	6822.0012
Sicherungen, Kit mit: Sicherung, 2 A, träge, 5 x 20 mm (10 St.) Sicherung, 4 A, träge, 5 x 20 mm (10 St.) Sicherung, 10 A, träge, 250V, 6,3 x 32 mm, 250 V (5 St.)	6820.0026
Signalkabel (6-pol. Mini-DIN)	6000.1004
Spritzen für WPS-3000(T)PL (RS) Nano/Cap und WPS-3000TBPL Nano/Cap 25μL 100 μL 250 μL, für analytische Anwendungen	6822.0001 6822.0002 6822.0003
Spritzen für WPS-3000TBPL Analytisch 100 μL 250 μL	6822.0002 6822.0003
Spritzen für WPS-3000TFC und WPS-3000TBFC 100 μL 250 μL Optimiert für Nano/Kap-Anwendungen	6822.0002 6822.0003
25μL	6822.0001

Bezeichnung	BestNr.
Spritzen-Bio-Kit (250 μL) (biokompatible Ausführung) für WPS-3000(T)PL (RS) Nano/Cap und WPS-3000TBPL Nano/Cap, enthält Spritze, 250 μL Nadel, 15 μL, Fused Silica, nanoViper Pufferschleife, 500 μL, PEEK/PTFE Probenschleife, 125 μL, PEEK	6821.0031
Spritzenkit (250 µL) für WPS-3000(T)PL (RS) Nano/Cap, enthält Spritze, 250 µL Nadel, 15 µL, Fused Silica, nanoViper Pufferschleife, 500 µL, SST/PTFE Probenschleife, 125 µL, nanoViper	6820.0031
Stator für Ventile mit 2-Positionen und 6 Ports Standardventil (Edelstahl) 900 bar-Ventil Biokompatibles Ventil (PAEK-Stator) Biokompatibles Ventil (Titan-Stator)* *Der Titan-Stator kann nicht für das Fraktionierventil verwendet werden.	6820.0012 6826.0012 6820.0034 6820.0049
Stator für Ventile mit 2-Positionen und 8 Ports (Micro Fraction Collection Option) Standardventil (Edelstahl) Biokompatibles Ventil	6820.0053 6821.0053
Transportfläschchen (zylindrisch, 10 mL), mit Kappe und Septum (5 Stück)	6820.0023
Transportfläschchenhalter (5 Positionen) für 10mL-Transportfläschchen	6820.4073
USB-Kabel, Typ A auf Typ B, High-Speed USB 2.0 (Kabellänge: 1 m)	6035.9035
USB-Kabel, Typ A auf Typ B, High-Speed USB 2.0 (Kabellänge: 5 m)	6911.0002
Vorstecher	6820.2400
Waschflasche WPS (2 Stück), einschließlich Halteklammer und Schlauch	6820.4075
Werkzeug zur Vormontage der biokompatiblen Fittinge	6000.0065

# 11 Index

Α

Anschluss	
Digital I/O	
Signalkabel	
Waschfllüssigkeit	
Anschlüsse	
Degaser	
elektrisch	35
Fluidik	27
Geräterückseite	22
Spritzenventil	55
USB	
Anzeige	21
Äquilibrieren	62
Auspacken	
Außerbetriebnahme	
Autosampler Drainage	
Autosamplerkonfigurationen	13
Autosampler-spezifische Informationen	

#### B

Back	77
Back-Flush-Modus	135, 138
Batch	73
Bedienelemente	21
Bedienungsanleitung	1
Beleuchtung	27
Betrieb	
Anzeige	21
Bedienelemente	21
Chromeleon	67
Einschalten	63
Empfehlungen	90
Sicherheitsmaßnahmen	4
Softkeys	74
User Interface	63
Biokompatible Autosampler	14, 146
- •	

#### С

Cancel	78
Chromeleon	29, 67
Autosampler installieren	40
Autosampler konfigurieren	41
Installationsprogramm	
Lizenz	29
Program Wizard	71
Server Configuration Program	38, 40
Server Monitor	38, 40
Service-Funktionen	91
Spezielle Funktionen	91
Steuerung	67
Steuerung automatisch	71
Steuerung direkt	68

USB-Treiber	
Chromeleon-Dialoge	
Error Levels	
General	
Inputs	
Options	
Relays	
Segments / Pump Link	
Sharing	
Commands (Dialogfenster)	
Configuration (Menü)	
Control (Menü)	79
Control Panel	
Coolant Temperature	
Cooling Power	
Cooling Pump Power	

## D

DCMSLink		48
Deep-Well-Plates		25
Degaser		52
Device View		72
Diagnostics (Menü)		81
Dialogfenster Commands		68
Digital I/O	23,	35
Dispense Destination		80
DispenseSpeed		85
Display		74
Display & Soft Keys		82
DispSpeed		85
Drainage		
Autosampler		56
System		56
DrawDelay		85
DrawSpeed		85
Drehteller		24
Konfiguration ändern		87
Tray Shake		94
Tray-Test		95
-		

## E

Einschalten	
Einstechtiefe (Vorstecher)	
Elektrischer Anschluss	
Entgasen	
Error Levels (Dialogfenster)	
Ersatzteile	
Execute	

#### F

Fan Power	
Fehlermeldungen	103, 104
Fehlersuche	
Fehlermeldungen	

LED blinkt	
Wellness (Chromeleon)	91
Firmware aktualisieren	
Firmware Download	
Fläschchenerkennung	
Fluidische Anschlüsse	
FlushVolume	
Fraction Collection	15
Fraktionierventil	
Fraktionierventil (WPS-3000FC)	
Fraktionssammlung	
Full-Loop-Injektion	
Funktionsprinzip	
WPS-3000FC	
WPS-3000PL	
Funktionstaste	
Back	
Cancel	
Clear	78.104
Execute	
Menu	
Next	78 104
Ok	78
Pfeil nach oben	77
Pfeil nach rechts	77
Pfeil nach unten	77
Prev	78 104
Prime	70, 101
Rotate	75
Select	
Wash	75
••••••••••••••••••••••••••••••••••••••	

#### G

Gefäßhöhen (Probengefäße)	25
General (Dialogfenster)	41
Gerätebeschreibung	
Biokompatible Autosampler	14, 146
Funktionsprinzip1	12, 135, 138
Konfigurationen	13
Temperaturregelung	15
Überblick	11
Gerätedisplay	
Helligkeit	
Kontrast	
Softkeys	74
Statusanzeige	64
Geräterückseite	
Digital I/O	23
Netzanschluss	23
Sicherung	23
USB	23
Geräterückseite (Anschlüsse)	
Netzkabel	
Signalkabel	
USB	

#### $\mathbf{H}$

Heat Sink Temperature
-----------------------

#### I

Inbetriebnahme	
Anschluss Waschflüssigkeit	
Äquilibrieren	
Kapillaren anschließen	
Inject Valve	80
Injektion (Synchronisation)	43. 86
Injektionsbefehl	
Injektionsmethode festelegen	
Injektionsmethoden	
Full-Loop	150
Low Dispersion	160
Mikroliter-Pickup	157
Partial-Loop	154
Injektionsparameter	
Injektionsventil	28, 125
WPS-3000FC	
WPS-3000PL	
Injizieren	
Innenansicht	
WPS-3000PL RS	17
WPS-3000TBFC	19
WPS-3000TBPL	17
WPS-3000TFC	19
Inputs (Chromeleon-Dialog)	46
Installation	
Anschlüsse (Geräterückseite)	
elektrischer Anschluss	
in DCMSLink	
Mikrotiterplatte	65
Mikrotiterplatten	58
Probenhalter	58, 65
Probenschleife	124
Pufferschleife	116
Reagensfläschchen	61
Spritzenwechsel	113
Standort	
Transportfläschchen	61
unter Chromeleon	
Verbinden	

## K

Kapillaren (Viper)	51
Kapillaren anschließen	51

## L

Leak Sensor Mode	
Leaksensor	
LoopOverfill	
Low Dispersion-Modus	

#### Μ

Main (Menü)	
Maintenance (Menü)	80
Menüs	76
Configuration	
Control	79
Diagnostics	
Main	
Maintenance	80
Micro Fraction Collection Option	145
Mikroliter-Pickup	157
Mikrotiterplatten	24
Einsetzen	
Empfehlung	25
Probenort	65
Mikrotiterplattenerkennung	65
Module Address	

#### Ν

Nadel waschen	
Nadelhöhe	93
Nadelpaar	118
Nadelpaar-Position	119
Netzschalter	23
Neue Hardware gefunden (Assistent)	
Nominal Sample Temperature	79
Normalphasen-Anwendungen (WPS-300	0FC)5
Normalphasen-Kit (WPS-3000TFC)	
-	

# 0

Ok	78
Online-Entgasung	52
Operational Qualification	95
Optionales Zubehör	
WPS-3000(T)PL	171
WPS-3000TBFC	178
WPS-3000TBPL	173
WPS-3000TBPL Analytisch	175
WPS-3000TFC	176
Options (Dialogfenster)	44

#### P

Panel Tabset	69
Partial-Loop-Injektion	154
Performance Qualification	95
Predictive Performance	91
Probenfläschchen	
Probengefäß	
Probenhalter	24
automatische Erkennung	95
Bestücken	
Einsetzen	
Nadelhöhe	93
Probenort	65
Tray Shake	94

Tray Test	95
Probennadel	
waschen	
Probenort	65
Probenposition	97
Probenraumbeleuchtung	
Probenschleife	16, 124
Probenthermostatisierung	
Probenvorbereitung	
Program Wizard	71
Programm	
Assistent	71
Benutzerdefiniert	
erstellen	
Herunterfahren	
Probenvorbereitung	
Shutdown	
Standby	
Programmstart	73
Pufferschleife	16, 116
Pump Link	
PunctureDepth	93

## R

Reagensfläschchen	
Einsetzen	61
Probenort	65
Real Temperature	81
Relative Probenposition	
Relays (Chromeleon-Dialog)	46
Reset to Factory Defaults	82
RMA-Nummer	111

## S

Sample Cooling	79
SampleHeight	
Segments / Pump Link (Dialogfenster)	
Select	77
Self Test	
Server Configuration Program	
Server Monitor	
Service	
Allgemein	111
Firmware aktualisieren	
Fraktionierventil	125
Injektionsventil	125
Leaksensor	112
Probennadel waschen	120
Probennadel wechseln	121
Probenschleife	124
Pufferschleife	116
Sicherheitsmaßnahmen	111
Sicherungswechsel	
Spritzenwechsel	113
Ventil Auseinanderbauen	127
Ventil Reinigen	126

Ventil Zusammenbauen	129
Vorstecher wechseln	123
Sharing (Dialogfenster)	42
Shutdown-Programm	99
Sicherheit	3
Sicherheitsmaßnahmen	4
Sicherungsschlitten	23
Sicherungswechsel	131
Signalkabel	35
Softkeys	74, 76
Spezifikation	163
Spritze	. 16, 113
Spritze Spülen	56
Spritzenventil	55
Spülen (Spritze)	56
Standby-Programm	99
Statusanzeige	64
Steuerung	
Automatisch	71
Chromeleon	67
Direkt	68
Symbole	3
Synchronisation (Injektion)	43, 86
Syringe	
System Wellness	
Systemdrainage	56

#### Т

Tastenfunktionen	
Technische Daten	
Temperaturregelung	
TransLiquidHeight	
Transportfläschchen	
Einsetzen	61
Probenort	65
TransVialPunctureDepth	
Tray	
Tray Shake	
Tray-Test	

#### U

Udp-Kommando	96
USB	
USB-Anschluss	
USB-Konfigurationsdatei	
User Interface	63

#### V

129

Verwendungszweck	8
Vials	25
Viper-Kapillaren	51
Volume	84
Vorbereitung	49
Allgemein	49
Anschluss Waschflüssigkeit	52
Äquilibrieren	62
Kapillaren anschließen	51
Reagensfläschchen	61
Spritze spülen	56
Spritzenventil	55
Transportfläschchen	61
Vorstecher	123
Vorstecher (Einstechtiefe)	93

#### W

Wartung	
Außerbetriebnahme	
Wartungsintervalle	
Wartungsintervalle	
Waschen (Probennadel)	
Waschflasche	
Waschflüssigkeit	
Wash	
WashSpeed	
WasteSpeed	85
Wechseln	
Probennadel	
Probenschleife	
Pufferschleife	
Sicherungen	
Spritze	
Vorstecher	
Wellness	
Well-Plates	
WPS-3000FC	
Betrieb	
Fraktionierventil	
Funktionsprinzip	
Injektionsventil	
Nano/Kap-Anwendungen	
Normalphasen-Anwendungen	5, 144
Normalphasen-Kit	
Service	144
WPS-3000PL	
Funktionsprinzip	
Micro Fraction Collection Option	
1	

# Z

Zubehör	
Optional	171
Standardzubehör	165