



UltiMate 3000 Series

Pump Series

Operating Instructions

(Original Instructions)



Revision: 1.3 Date: November 2011

© 2011 Thermo Fisher Scientific Doc No. 4820.3551

CE

Declaration of Conformity

(Original Declaration)

Product: UltiMate 3000 Series Pump

Types: LPG-3400AB, LPG-3400M and LPG-3400MB DGP-3600A, DGP-3600AB, DGP-3600M and DGP-3600MB HPG-3200P

Dionex Softron GmbH herewith declares conformity of the above products with the respective requirements of the following regulations:

- Machinery Directive 2006/42/EC
- EMC Directive 2004/108/EC

The safety of the machinery was evaluated based on the following standard:

 EN ISO 12100:2010
 Safety of machinery - General principles for design Risk assessment and risk reduction

The electrical safety of the products was evaluated based on the following standard:

• DIN EN 61010-1:2002 Safety requirements for electrical equipment for measurement, control and laboratory use, Part 1: General Requirements

The Electromagnetic Compatibility (EMC) of the products was evaluated based on the following standard:

• DIN EN 61326: 2006 Electrical equipment for measurement, control and laboratory use EMC Requirements

The protection requirements specified in the low-voltage directive 2006/95/EC are met.

Responsible for the technical CE documentation is the manufacturer (see further down).

This declaration is issued for the manufacturer

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

by the President, Dr. Peter Jochum. November 2, 2011

| Certificate no. | CU 72051793 02 | |
|--|--|---|
| License Holder: Dionex Softro Dornierstr. 4 | n GmbH Dionex Sof Dornierstr | Plant: tron GmbH . 4 |
| 82110 Germeri Germany | ng 82110 Germ Germany | mering |
| Test report no.: U | SA-RC 30581482 002 Client Reference | e: Burkhard Seyferth |
| Tested to: U | L 61010-1:2004 AN/CSA-C22.2 61010-1:2004 | |
| Certified Product: | HPLC Pump | License Fee - Units |
| Certified Product: Addition: Model Designa | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z | License Fee - Units = B, blank) |
| Certified Product: Addition: Model Designa Appendix: 1, | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 | License Fee - Units = B, blank) |
| Certified Product: Addition: Model Designa Appendix: 1, | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 | License Fee - Units = B, blank) |
| Certified Product: Addition: Model Designa Appendix: 1, | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 | License Fee - Units = B, blank) |
| Certified Product: Addition: Model Designa Appendix: 1, | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 | License Fee - Units = B, blank) |
| Certified Product: Addition: Model Designa Appendix: 1, | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 | License Fee - Units = B, blank) |
| Certified Product: Addition: Model Designa Appendix: 1, Licensed Test mark | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 Sig | License Fee - Units = B, blank) natures Date of Issue (day/mo/yr) 20/12/2006 |
| Certified Product: Addition: Model Designa Appendix: 1, Licensed Test mark | HPLC Pump ation: ISO-3100yz, LPG-3400yz, DPG-3600yz, HPG-3X00yz (X = 2, 4; y = M, A, P; z 1-2 1-2 Sign J. Sign Stephan Schmitt | License Fee - Units = B, blank) natures Date of Issue (day/mo/yr) 20/12/2006 |

Table of Contents

| 1 | In | ntroduction | 1 |
|---|-----------|--|----|
| | 1.1 | How to Use This Manual | 1 |
| | 1.2 | Safety Information | 2 |
| | 1.2 | 2.1 Symbols on the Instrument and in the Manual | 2 |
| | 1.2 | 2.2 General Safety Precautions | |
| | 1.2 | 2.3 Consignes Générales de Sécurité | 5 |
| | 1.3 | Intended Use | 8 |
| | 1.4 | Federal Communications Commission (FCC) Note | |
| 2 | O | verview | 9 |
| _ | 21 | Unit Description | 0 |
| | 2.1 | Operating Principle | |
| | 2.2 | Pump Configurations | 10 |
| | 2.5 ?? | 3 1 Overview | 12 |
| | 2.2 | 3 2 Biocompatible Pumps | 12 |
| | 2.4 | Interior Components | 16 |
| | 2.5 | Front Panel Flements | 10 |
| | 2.5 | Rear Panel | 17 |
| | 2.0 | 6.1 Power Switch | |
| | 2.6 | 6.2 Fuse Cartridge | |
| | 2.6 | 6.3 USB Connector | 19 |
| | 2.6 | 6.4 Pressure (Analog Pressure Output) | |
| | 2.6 | 6.5 Digital I/O | |
| | 2.6 | 6.6 Solvent Rack | |
| | 2.7 | Fluid Connections | |
| | 2.8 | Rear Seal Wash System | |
| | 2.9 | Outlet Unit | |
| | 2.9 | 9.1 Inline Filter | |
| | 2.9 | 9.2 Mixing Chamber | |
| | 2.10 | Leak Sensor | |
| | 2.11 | Vacuum Degasser | |
| | 2.12 | Chromeleon Software | |
| | 2.13 | System Wellness, Predictive Performance, and Diagnostics | |
| 3 | In | nstallation | |
| | 3.1 | Facility Requirements | |
| | 3.2 | Unpacking | |
| | 33 | Positioning the Pump in the UltiMate 3000 System | 30 |
| | 3.4 | Connecting the Pump | 32 |
| | 3.4 | 4.1 General Information | |
| | 3.4 | 4.2 Connecting the USB Cable | |

| 3.4.3 | Connecting the Power Cord | |
|--|--|---|
| 3.4.4 | Connecting the Solvent Rack and Digital I/O | |
| 3.5 Se | ting Up the Pump in Chromeleon | |
| 3.5.1 | Loading the USB Driver for the Pump | |
| 3.5.2 | Installing the Pump | |
| 3.5.3 | Configuring the Pump | |
| 3.5 | .3.1 Initial Installation | |
| 3.5 | .3.2 Changing the Configuration Properties | 44 |
| 4 Prepa | ration for Operation (Startup) | 45 |
| 4.1 Ov | erview of Actions | 45 |
| 4.2 Ge | neral Precautions for Connecting Capillaries | |
| 4.3 So | lvent Reservoirs | 47 |
| 4.3.1 | General Precautions | 47 |
| 4.3.2 | Connecting the Solvent Reservoirs | |
| 4.4 Co | nnecting the Rear Seal Wash System | 49 |
| 4.4.1 | Filling the Internal Wash Liquid Reservoir | 50 |
| 4.4.2 | Installing and Filling the External Wash Liquid Reservoir | |
| 4.5 Co | nnecting the Drain Tubing | 53 |
| 4.6 Pu | rging the Pump | 54 |
| 4.6.1 | Purging the Pump Manually | 54 |
| 4.6.2 | Purging the Pump via the WPS-3000SL Autosampler | |
| | | |
| 4.7 Eq | uilibrating the System | 57 |
| 4.7 Eq 5 Oper | ation and Maintenance | 57 |
| 4.7 Eq 5 Oper | ation and Maintenance | 57 59 |
| 4.7 Eq 5 Oper 5.1 Po | ation and Maintenance wer-Up | 57 59 59 |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta | ation and Maintenance | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op | ation and Maintenance wer-Up tus Screens eration with Chromeleon | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3 2 | ation and Maintenance | 57 59 60 61 61 61 |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 | ation and Maintenance | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di | ation and Maintenance wer-Up tus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4 1 | ation and Maintenance wer-Up itus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control Splay Screens (Function Keys and Menus) Showing the Function Keys | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 | ation and Maintenance wer-Up tus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control Splay Screens (Function Keys and Menus) Showing the Function Keys Pump Menus | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 | ation and Maintenance wer-Up itus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys Pump Menus .2.1 General Menu Layout and Structure | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up itus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys Pump Menus .2.1 General Menu Layout and Structure. .2.2 Main Menu | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up tus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys Pump Menus .2.1 General Menu Layout and Structure .2.2 Main Menu .2.3 Control Menu | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up tus Screens eration with Chromeleon connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys Pump Menus .2.1 General Menu Layout and Structure .2.2 Main Menu .2.3 Control Menu .2.4 Preferences Menu | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up itus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys. Pump Menus .2.1 General Menu Layout and Structure .2.2 Main Menu .2.3 Control Menu .2.4 Preferences Menu .2.5 Diagnostics Menu | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up tus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys. Pump Menus .2.1 General Menu Layout and Structure. .2.2 Main Menu .2.3 Control Menu .2.4 Preferences Menu .2.5 Diagnostics Menu .2.6 Configuration Menu | 57 59 60 61 62 65 67 68 69 70 71 72 73 74 |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up tus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control splay Screens (Function Keys and Menus) Showing the Function Keys Pump Menus .2.1 General Menu Layout and Structure .2.2 Main Menu .2.3 Control Menu .2.4 Preferences Menu .2.5 Diagnostics Menu .2.6 Configuration Menu erational Settings Choosing the Solvents | |
| 4.7 Eq 5 Oper 5.1 Po 5.2 Sta 5.3 Op 5.3.1 5.3.2 5.3.3 5.4 Di 5.4.1 5.4.2 5.4 5.4 5.4 5.4 5.4 5.4 5.4 5.4 | ation and Maintenance wer-Up tus Screens eration with Chromeleon Connecting to Chromeleon Direct Control Automated Control Splay Screens (Function Keys and Menus) Showing the Function Keys. Pump Menus .2.1 General Menu Layout and Structure. .2.2 Main Menu .2.3 Control Menu .2.4 Preferences Menu .2.5 Diagnostics Menu .2.6 Configuration Menu erational Settings Choosing the Solvents Linking the Pump to the Autosampler | |

| | 5.5.4 | Setting the Flow Rate, Flow Acceleration, and Flow Deceleration | 78 |
|-----|--------------------|--|-----|
| | 5.5.5 | Setting the Pressure Limits | 79 |
| | 5.5.6 | Recording the Pump Pressure | 80 |
| | 5.5.7 | Activating Rear Seal Washing | 81 |
| | 5.5.8 | Purging the Pump | |
| | 5.5.9 | Activating Leak Detection | 84 |
| | 5.5.10 | Adjusting the Screen Brightness or Contrast | 85 |
| | 5.5.11 | SmartStartup and SmartShutdown | 85 |
| 5.0 | 6 Sp | ecial Chromeleon Functions | 86 |
| | 5.6.1 | Predictive Performance | 86 |
| | 5.6.2 | Pump Diagnostics | 87 |
| | 5.6.3 | Setting a Gradient Curve | 88 |
| | 5.6.4 | Using the Digital Inputs and Outputs (Digital I/O) | 88 |
| | 5.6.5 | Operational Qualification und Performance Qualification | 89 |
| | 5.6.6 | Enabling the Double Flow Mode | 89 |
| 5.′ | 7 Sh | atting Down the Pump | 90 |
| 5.8 | 8 Ro | utine and Preventive Maintenance Intervals | 92 |
| | T I | | 0.5 |
| 6 | Trou | oleshooting | 95 |
| 6. | 1 Ov | erview | 95 |
| 6.2 | 2 Me | ssages on the Front Panel Display | 96 |
| 6. | 3 Ch | romeleon Diagnostics Messages | 101 |
| 6.4 | 4 Op | erating Problems | 103 |
| | 1 | | |
| 7 | Servi | ee | 109 |
| 7. | 1 Ge | neral Notes | 109 |
| 7.2 | 2 Dr | ving the Leak Sensor | 110 |
| 7 | 3 Fil [.] | er Holder and Filter Frit in the Inline Filter | 111 |
| / | 731 | Checking the Permeability of the Filter Frit | 112 |
| | 7.3.2 | Exchanging the Filter Holder or Filter Frit. | 112 |
| 74 | 4 Re | placing the Check Valves | 113 |
| 7.4 | 5 Pis | tons and Piston Seals | 114 |
| / | 751 | Visually Inspecting the Pump for Piston Seal Leakage | 117 |
| | 752 | Replacing the Piston Seals | 119 |
| | 7.5 | 2.1 Removing the Pump Head and Piston | 120 |
| | 7.5 | 2.2 Removing the Piston Seals and Support Ring | 123 |
| | 7.5 | .2.3 Cleaning the Pistons | 124 |
| | 7.5 | .2.4 Installing the Piston, Piston Seals, Support Ring, and Pump Head | 125 |
| | 7.5 | .2.5 Recommended Actions after Exchanging the Piston Seals | 130 |
| 7.0 | 6 Re | placing the Fuses | 131 |
| 7. | 7 Va | cuum Degasser | 132 |
| 7 9 | 8 Te | ting the Pump for Leakage | 133 |
| /.(| | and the second for th | 155 |

| 8 O | ptimizing the Pump for Special Application | |
|------|---|------------|
| 8.1 | Mixing Chamber Extension | |
| 8.2 | Micro Flow Kit | |
| 8.3 | Static Mixer | 141 |
| 9 Ir | nside Front Panel Views and Fluid Connections | 143 |
| 9.1 | Isocratic Pump (IS0-3100A) | |
| 9. | 1.1 Inside Front Panel | |
| 9. | 1.2 Fluid Connections | |
| 9.2 | Low-Pressure Gradient Pumps | |
| 9. | 2.1 LPG-3400 | |
| | 9.2.1.1 Inside Front Panel | 146 |
| | 9.2.1.2 Fluid Connections | 147 |
| 9. | 2.2 DGP-3600 | |
| | 9.2.2.1 Inside Front Panel | |
| | 9.2.2.2 Fluid Connections | 149 |
| 9.3 | High-Pressure Gradient Pumps | |
| 9. | 3.1 HPG-3200 | |
| | 9.3.1.1 Inside Front Panel | |
| 0 | 9.3.1.2 Fluid Connections | |
| 9. | 3.2 HPG-3400 | |
| | 9.3.2.1 Inside Front Panel | |
| 0 | 9.5.2.2 Fluid Collifections | 134 156 |
| 9. | 9.3.3 Supported Gradient Combinations | 150 156 |
| | 9.3.3.2 Double Flow Mode | 150 |
| 9. | 3.4 HPG-3200P Semipreparative Pump | |
| 10 T | echnical Information | |
| 11 A | ccessories, Spare Parts, and Consumables | |
| 11.1 | Standard Accessories | |
| 11.2 | 2 Optional Accessories | |
| 11.3 | Consumables and Spare Parts | |
| 12 R | eference Information | |
| 12.1 | Chemical Resistance of PEEK | 175 |
| 12.1 | Solvent Miscibility | 178 |
| 12.2 | Properties of Common Solvents | 170 |
| 12.3 | Safaty Information about Elammable Salvanta | 1/9 100 |
| 12.4 | sancey mormation about Frammable Solvents | |

| 13 App | endix | |
|---------|--|--|
| 13.1 | Connections in the Low- and High-Pressure Sections | |
| 13.2 | Manual Injection Valve | |
| 13.3 | Pin Assignment | |
| 13.3 | 1 D-Sub I/O (Digital I/O) | |
| 13.3 | 2 Solvent Rack | |
| 14 Inde | X | |

1 Introduction

1.1 How to Use This Manual

The layout of this manual is designed to provide quick reference to the sections of interest to the reader. However, in order to obtain a full understanding of the pump, Thermo Fisher Scientific recommends that you review the manual thoroughly before beginning operation.

Almost all descriptions in the manual apply to all pump types in the UltiMate[™] 3000 series pumps and cover both the standard (stainless steel) and biocompatible pumps. Therefore, the term "the pump" is used throughout the manual. If some detail applies to only one type or model, the model is identified by name. If only the pump name, for example DGP-3600, is used, the information applies to all pump models (DGP-3600A, DGP-3600M, and DGP-3600MB).

Notes: The device configuration may vary; therefore, not all descriptions necessarily apply to your particular instrument.

The descriptions in this manual refer to firmware version 3.05 and Chromeleon[™] 6.80 Service Release 11. If you want to operate the pump with Chromeleon 7, note the information on page 26.

This manual is provided "as is." Every effort has been made to supply complete and accurate information and all technical specifications have been developed with the utmost care. The information contained in this manual should not be construed as a commitment by Thermo Fisher Scientific. Thermo Fisher Scientific assumes no responsibility for any errors that may appear in this document. This document is believed to be complete and accurate at the time of publication. In no event shall Thermo Fisher Scientific be liable for incidental or consequential damages in connection with or arising from the use of this document. The information contained in this document is subject to change without notice.

All rights reserved, including those for photomechanical reproduction and storage on electronic media. No part of this publication may be copied or distributed, transmitted, transcribed, stored in a retrieval system, or transmitted into any human or computer language, in any form or by any means, electronic, mechanical, magnetic, manual, or otherwise, or disclosed to third parties without the express written permission of Thermo Fisher Scientific Inc.

Trademarks

Analyst is a registered trademark of AB Sciex. Compass and Hystar are trademarks of Bruker Daltronics. Empower is a trademark of Waters Corp. PEEK is a trademark of Victrex PLC. PharMed is a registered trademark of Saint -Gobain Performance Plastics. Windows and Windows Vista are registered trademarks of Microsoft Corp. All other trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries.

1.2 Safety Information

The CE Mark label and cTUVus Mark safety label on the instrument indicate that the instrument is compliant with the related standards (\rightarrow pages I and I).

1.2.1 Symbols on the Instrument and in the Manual

The table shows the symbols used on the instrument:

| Symbol | Description |
|--------------|---|
| ~ | Alternating current—Courant alternatif |
| - 0 | Power supply is on $(-)$ — L'instrument est mis sous tension $(-)$ and Power supply is off (\mathbf{O}) — L'instrument est mis hors tension (\mathbf{O}) |
| \triangle | Refer to the <i>Operating Instructions</i> to prevent risk of harm to the operator and to protect the instrument against damage. Référez-vous à ce manuel pour éviter tout risque de blessure à l'opérateur et/ou protéger l'instrument contre tout dommage. |
| 1 (1) | Label according to the "Measures for Administration of the Pollution Control of Electronic Information Products" (China RoHS) guideline Étiquette "Measures for Administration of the Pollution Control of Electronic Information Products" (China RoHS) |
| | WEEE (Waste Electrical and Electronic Equipment) label—For more information, see the WEEE Information section in the "Installation and Qualification Documents for Chromatography Instruments" binder. Étiquette WEEE (Waste Electrical and Electronic Equipment) —Pour plus d'informations, référez-vous au chapitre WEEE Information dans le classeur "Installation and Qualification Documents for Chromatography Instruments". |

At various points throughout the manual, messages of particular importance are indicated by certain symbols:

| i | Tip: | Indicates general information, as well as information intended to optimize the performance of the instrument. |
|------|----------------|--|
| Ŵ | Important: | Indicates that failure to take note of the accompanying information could cause wrong results or may result in damage to the instrument. |
| Ŵ | Important: | Indique que ne pas tenir compte de l'information jointe peut conduire à de faux résultat ou endommager l'instrument. |
| STOP | Warning: | Indicates that failure to take note of the accompanying information may result in personal injury. |
| STOP | Avertissement: | Indique que ne pas tenir compte de l'information jointe peut entraîner des blessures corporelles. |

1.2.2 General Safety Precautions

When working with analytical instrumentation, you should know the potential hazards of using chemical solvents. Wear appropriate protective clothing.

I Tips: Before initial operation of the pump, make sure that you are familiar with the contents of this manual.

Observe any warning labels on the device and see the related sections in these *Operating Instructions*.

For the general safety precautions in French, see page 5.

To avoid the possibility of personal injury and damage to the instrument, observe the following general safety precautions when operating the instrument or carrying out any maintenance work:

- Install the HPLC system in a well-ventilated laboratory. If the mobile phase includes volatile or flammable solvents, do not allow them to enter the workspace.
- For minimum interference effects, all components of the analytical system should be connected to the same mains output (same phase).
- The pump is primed with 2-propanol. During initial operation of the pump, make sure that the solvents used are miscible with 2-propanol. Otherwise, follow the appropriate intermediate steps.
- The front panel tilts upward. To prevent damage to the pump when lifting or moving, always lift by the bottom or sides of the unit.
- Do not place any heavy objects on the open front panel door. This may damage the door.
- Always set a lower pressure limit for the HPLC pump. This prevents damage resulting from leakage or from running the pump dry.
- To avoid that the pressure calibration of the pump is impaired, turn on the pump only when the pump pressure is down. To ensure that the pressure is down, open the purge screw before turning on the pump.
- Never run the pump dry. Damage to the pistons or the piston seals could result.
- Thermo Fisher Scientific advises against recycling the solvents. This may impair the performance of the seals.
- When connecting the capillaries, make sure that the connectors are free from contaminants. Even minute particles may cause damage to the system (for example, flow splitter, flow control valve, and column).
- After operation, rinse out buffers and solutions that form peroxides.

- Before switching from buffer to organic solution, rinse the pump thoroughly with deionized water.
- When switching to another solvent, ensure that the new solvent is miscible with the one contained in the pump. Otherwise, the pump can be damaged; for example, by flocculation!
- If the pump flow is interrupted for longer periods (> 1 hour), turn off the lamps in any UV or RF detector connected to the pump. This will prevent evaporation in the flow cell.
- If you use solvents with a high salt content, do not operate the pump without rear seal washing (→ page 49) for a longer time (> 5 minutes). This may cause damage to the piston seals and the piston. Regularly exchange the liquid in the liquid reservoir of the rear seal wash system (at least once a week).
- Always use the frits recommended by Thermo Fisher Scientific to prevent particulate matters from entering the HPLC system. Using other frits may considerably affect the system performance.
- Do not use stainless steel frits with biocompatible versions of the pump. This renders he biocompatibility void.
- If the mobile phase includes volatile or flammable solvents, avoid open flames and sparks.
- If a leak occurs, turn off the instrument and remedy the situation immediately.
- When the panels are removed, dangerous electrical connections will be exposed. Disconnect the pump from all power sources before removing the panels. The enclosure should be opened by authorized service personnel only.
- Always replace blown fuses with original Dionex spare part fuses (\rightarrow page 131).
- Replace faulty power cords and communication cables.
- Many organic solvents and buffers are toxic. Know the toxicological properties of all mobile phases that you are using.
- The toxicological properties of many samples may not be well known. If you have any doubt about a sample, treat it as if it contains a potentially harmful substance.
- Wear goggles when handling mobile phases or operating the instrument. An eyewash facility and a sink should be close to the unit. If any mobile phase splashes on the eyes or skin, wash the affected area and seek medical attention.
- Dispose of waste mobile phase in an environmentally safe manner that is consistent with all local regulations. Do not allow flammable or toxic solvents to accumulate. Follow a regulated, approved waste disposal program. Never dispose of flammable or toxic solvents through the municipal sewage system

- Use only standard solvents (HPLC grade) and buffers that are compatible with all parts that may be exposed to solvents (\rightarrow page 159).
- In an UltiMate 3000 system, some components are made of PEEK[™]. While this polymer has superb chemical resistance to most organic solvents, it tends to swell when in contact with trichlormethane (CHCl₃), dimethyl sulfoxide (DMSO), or tetrahydrofuran (THF). In addition, it is attacked by concentrated acids, such as, sulfuric acid and nitric acid or a mixture of hexane, ethyl actate, and methanol. (Swelling or attack by concentrated acids is not a problem with brief flushing procedures.)
- Do not use PEEK tubing that is stressed, bent, or kinked.
- Before interrupting operation for several days or more or when preparing the pump for transport, observe the precautions for shutting down the pump (\rightarrow page 90).
- Use original Dionex spare parts only. Substituting non-Dionex parts or using non-Dionex accessories may impair the performance of the instrument.
- Do not use the pump in ways other than those described in this manual.

1.2.3 Consignes Générales de Sécurité

Veuillez noter: Avant de commencer à utiliser la pompe, assurez-vous que vous vous êtes familiarisés avec le contenu de ce manuel.

Observez des étiquettes d'avertissement sur l'appareil et référez-vous aux sections correspondantes dans ce mode d'emploi.

Veuillez observer les consignes générales de sécurité suivantes lorsque vous utilisez l'instrument ou que vous procédez à des opérations de maintenance.

- Installez le système HPLC dans un laboratoire bien ventilé. Si la phase mobile contient des solvants volatils ou inflammables, empêchez qu'ils ne pénètrent dans l'espace de travail.
- Afin d'éviter au maximum les interférences, tous les éléments du système analytique doivent être raccordés à la même ligne secteur (même phase).
- La pompe est stockée sous 2-propanol. Au cours démarrage de la pompe, assurez-vous que les solvants utilisés soient miscibles avec le 2-propanol. Sinon, suivez les étapes intermédiaires appropriées.
- Le panneau avant bascule vers le haut. Afin d'éviter d'endommager la pompe lorsque que vous la soulevez ou la déplacez, saisissez-la toujours par le bas ou les côtés de l'unité.
- Ne placez aucun objet lourd sur la porte ouverte du panneau avant. Ceci pourrait endommager la porte.

- Réglez toujours une limite de pression minimum pour la pompe HPLC. Ceci prévient les dommages résultant de fuites ou du fonctionnement à sec de la pompe.
- Afin d'éviter que le calibrage de pression de la pompe ne soit pas entravé, mettez en marche la pompe seulement quand le pompe est sans pression. Toujours ouvrez la vis de purge avant mettre la pompe en marche.
- Ne faites jamais fonctionner la pompe à sec. Il peut en résulter des dommages aux pistons ou aux joints de piston.
- Thermo Fisher Scientific déconseille de recycler les solvants. Ceci peut nuire aux performances des joints.
- Lorsque vous connectez les capillaires, assurez-vous que les raccords sont exempts de tout contaminant. Même d'infimes particules peuvent causer des dommages au système (ex. diviseur de débit, vanne de régulation de débit et colonne).
- Après utilisation, purgez le système des tampons et des susceptibles de former des peroxydes.
- Lorsque vous passez d'une solution saline à un solvant organique, effectuez un rinçage intermédiaire de la pompe à l'eau dé-ionisée.
- Lorsque vous passez à un autre solvant, assurez-vous que le nouveau solvant soit miscible avec celui qui se trouve dans la pompe. Dans le cas contraire, la pompe peut être endommagée; par exemple, par des floculations!
- Si le débit de la pompe est interrompu pour des périodes prolongées (> 1 heure), éteignez les lampes de tout détecteur UV ou RF raccordé à la pompe. Ceci empêchera l'évaporation dans la cellule.
- Si vous utilisez des phases mobiles avec une forte teneur en sel, ne faites pas fonctionner la pompe sans rinçage du joint arrière pendant un temps prolongé (> 5 minutes). Ceci peut endommager les joints de piston et le piston (→ page 49). Remplacer régulièrement le liquide dans le réservoir du système de rinçage du joint arrière (au moins une fois par semaine).
- Utilisez toujours les frittés recommandés par Thermo Fisher Scientific afin d'empêcher les particules étrangères d'entrer dans le système HPLC. Utiliser d'autres frittés peut affecter considérablement les performances du système.
- N'employez pas des frittes d'acier inoxydable avec des versions biocompatible de la pompe. Ceci rend vide de compatibilité biologique de la pompe.
- Si la phase mobile contient des solvants volatils ou inflammables, évitez les flammes nues et les sources d'étincelles à proximité.
- Si une fuite survient, arrêtez l'instrument et résolvez le problème immédiatement.

- Quand les capots de protection de l'appareil sont démontés, vous êtes exposés à des connexions électriques sous haute tension deviennent accessibles. Débranchez le passeur d'échantillon de toute source d'alimentation électrique avant de retirer les capots. Ne démontez les capots de protection que si cela est explicitement demandé au cours de ces instructions.
- Remplacez toujours les fusibles grillés par des fusibles de rechange d'origine Dionex (→ page 131).
- Remplacez les cordons d'alimentation électrique et les câbles de communication défectueux.
- De nombreux solvants organiques et solutions salines sont toxiques. Informez-vous des propriétés toxicologiques de toutes les phases mobiles que vous utilisez. Portez le vêtement de protection approprié.
- Les propriétés toxicologiques de nombreux échantillons peuvent être mal connues. Au moindre doute concernant un échantillon, traitez-le comme s'il contenait une substance potentiellement dangereuse.
- Portez des lunettes de protection lorsque vous manipulez des phases mobiles ou que vous utilisez l'instrument. Une installation permettant de se laver les yeux ainsi qu'un lavabo doivent se trouver à proximité du système. Si une phase mobile, quelle qu'elle soit, entre en contact avec vos yeux ou votre peau, rincez abondamment la zone affectée à l'eau, puis.
- Débarrassez-vous de tous les déchets de phase mobile de manière écologique, conformément à la règlementation en vigueur au niveau local. Empêchez impérativement l'accumulation de solvants inflammables et/ou toxiques. Suivez un programme d'élimination des déchets règlementé et approuvé. Ne jetez jamais de solvants inflammables et/ou toxiques dans le système municipal d'évacuation des eaux usées.
- Utilisez uniquement des solvants (qualité HPLC) et des solutions salines compatibles avec les matériaux exposés phase mobiles (→ page 159).
- Dans un système UltiMate 3000, certaines composantes sont en PEEK. Bien que ce polymère présente une excellente résistance chimique à la plupart des solvants organiques, il a tendance à gonfler lorsqu'il est en contact prolongé avec du chloroforme (CHCl₃), du diméthyle sulfoxyde (DMSO) ou du tétrahydrofurane (THF). De plus, il est attaqué par des acides concentrés tels que l'acide sulfurique et l'acide nitrique ou d'un composé du hexane, éthyle acétate et méthanol. (Ces acides peuvent cependant être utilisés dans le cadre de procédures de nettoyage, à condition que l'exposition soit brève.)
- N'utilisez pas de tubes PEEK écrasés, pliés ou abimés.
- Avant d'interrompre le fonctionnement pendant plusieurs jours ou plus, observez les précautions figurant en Shutting Down the Pump (→ page 90).

- Utilisez des pièces de rechange d'origine Dionex. Effectuer des remplacements par des pièces ne provenant pas de Thermo Fisher Scientific ou utiliser des accessoires ne provenant pas de Thermo Fisher Scientific peut affecter les performances de l'instrument.
- N'utilisez pas la pompe de manière autre que celles décrites dans ce manuel.

1.3 Intended Use

The pump is designed to perform equally well as a dependable system for routine analyses or as a sophisticated research instrument for use in HPLC (high performance liquid chromatography) applications, especially as part of the UltiMate 3000 system. However, it can also be used with other HPLC systems if adequate control inputs and outputs are available. A PC with USB port is required.

The pump is operated with the Chromeleon Chromatography Management System. Being part of the UltiMate 3000 system, the pump can also be operated with other data systems, such as

• Analyst[®], CompassTM/HyStarTM or XcaliburTM

To do so, installation of the DCMSLink (Dionex Chromatography Mass Spectrometry Link) software is required in addition to the installation of the data system.

 EmpowerTM Installation of the Dionex Instrument Integration Software is required in addition to the installation of the data system.

For more information, contact the Thermo Fisher Scientific sales organization for Dionex HPLC Products.

Please note that the pump may be operated only using the accessories originally supplied with the units (\rightarrow page 163) and within their technical specifications (\rightarrow page 158).

If there is any question regarding appropriate usage, contact Thermo Fisher Scientific before proceeding.

Thermo Fisher Scientific cannot be held liable for any damage, material or otherwise, resulting from inappropriate or improper use of the instrument.

1.4 Federal Communications Commission (FCC) Note

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the U.S. FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference, in which case the user will be required to correct the interference at his expense.

2 Overview

2.1 Unit Description

The pump is a modern high-quality instrument designed for HPLC analysis, especially as part of the UltiMate 3000 system. The pump was especially developed for routine HPLC analysis and can be used in numerous laboratory environments. The instrument performs equally well as a flexible and reliable module for routine analysis and sophisticated research tasks:

- The patented isokinetic pre-compression allows a precise and almost pulse-free flow.
- The technical specification meets the highest requirements for flow rate reproducibility, zero pulsation, and operational reliability.
- All pumps (except the semipreparative pump) are fitted with floating pistons, allowing compensation for small mechanical tolerances within the specification and thus enhancing the robustness of the pump.
- Various monitoring and diagnostic features are provided for optimum system performance and reliability (\rightarrow page 27).
- For the secure and functional positioning of the solvent reservoirs on top of the pump, the Solvent Racks of the UltiMate 3000 series are available from Thermo Fisher Scientific (→ page 14). Except for the SR-3000, all Solvent Racks include an integrated vacuum degasser.
- The pump can be fully controlled by the Chromeleon Chromatography Management System, providing a high degree of system integration.
- All parts that may be exposed to solvents are made of materials that provide optimum resistance to the most commonly used solvents and buffer solutions.
- The UltiMate 3000 pump series includes also biocompatible pump versions. For information about the characteristics of these pumps, see page 14.

2.2 Operating Principle

The pump is a zero-pulsation, serial dual-piston pump with electronic compressibility compensation. The two pump heads are connected in series. The solvent passes through both pump heads—working and equilibration head—successively.

Continuous delivery is achieved as follows: The working head delivers at the appropriate flow rate while simultaneously filling the serially connected equilibration head. The latter serves as a reservoir and delivers while the working head carries out the suction stroke.

The characteristic feature of the patented isokinetic pre-compression is the 120-degree overlapping phase of the delivery strokes of the working and equilibration heads. When delivering compressible liquids without controlled pre-compression, the pulsation increases as the operating pressure increases, since part of the delivery stroke is required for compressing the solvent in the pump head.

Pulsation during the pre-compression phase is reduced to a minimum by velocity modulation of the drive. The highly constant delivery is ensured by a patented secondary control system (automatic compressibility compensation). The flow rate is always kept constant in relation to the atmospheric pressure.

The pictures illustrate how the pump operates:



Fig. 1: LPG-3400 pump with integrated vacuum degasser



Fig. 2: DGP-3600 pump



Fig. 3: HPG-3200 pump

For information about the pump configurations and options, see page 12.

2.3 Pump Configurations

2.3.1 Overview

A binary high-pressure gradient system, optionally equipped with the "2 from 4" solvent selector, is available in addition to an isocratic analytical pump and the quaternary low-pressure gradient pump with integrated degasser. For semipreparative applications, a binary high-pressure gradient pump with a maximum flow of 100 mL/min is provided.

A dual low-pressure gradient pump completes the pump series. You can use this pump to run two ternary gradients independently from each other. The dual gradient pump incorporates two independent pumps in a single housing. Thus, you can use two pumps without occupying additional bench space.

A dynamic mixing chamber with variable volume that can be adjusted to individual requirements offers additional flexibility.

| Pump | Description | Part No. | Option |
|--|--|-----------|---|
| ISO-3100A | Isocratic pump (analytical; 1 solvent) | 5035.0010 | |
| LPG-3400A Low-pressure gradient pump (analytical; 4 solvents) with integrated vacuum degasser and mixing chamber | | 5035.0015 | Mixing Chamber Extension Kit (→ page 136) Micro Flow Kit (→ page 138) |
| LPG-3400M | Low-pressure gradient pump optimized for micro flows (4 solvents) with integrated vacuum degasser. The pump has no mixing chamber. | 5035.0045 | |
| LPG-3400AB | Same as LPG-3400A, but biocompatible version | 5037.3015 | Mixing Chamber Extension Kit (→ page 136) Micro Flow Kit (→ page 138) |
| LPG-3400MB | Same as LPG-3400M, but biocompatible version | 5037.0055 | |
| DGP-3600A | Dual low-pressure gradient pump (analytical): Two separate pumps with integrated mixing chambers in one enclosure (2x3 solvents) | 5035.0014 | Mixing Chamber Extension Kit (→ page 136) Micro Flow Kit (→ page 138) |
| DGP-3600M | Dual low-pressure gradient pump optimized for micro flows: Two separate pumps are installed in one enclosure (2x3 solvents). The pumps have no mixing chambers. | 5035.0050 | |

The pump is available in the following configurations:

| Pump | Description | Part No. | Option |
|------------|---|-----------|---|
| DGP-3600AB | Same as DGP-3600A, but biocompatible version | 5037.0014 | Mixing Chamber Extension Kit (→ page 136) Micro Flow Kit (→ page 138) |
| DGP-3600MB | Same as DGP-3600M, but biocompatible version | 5037.0060 | |
| HPG-3200A | High-pressure gradient pump (analytical; 2 solvents) with integrated mixing chamber | 5035.0016 | Mixing Chamber Extension Kit $(\rightarrow page 136)$ |
| HPG-3200M | High-pressure gradient pump optimized for micro flows (2 solvents). The pump has no mixing chamber. | 5035.0018 | |
| HPG-3200P | High-pressure gradient pump (semipreparative; 2 solvents) with integrated mixing chamber and mixing chamber extension | 5035.0025 | Mixing Chamber Extension Kit (→ page 136) |
| HPG-3400A | High-pressure gradient pump (analytical) with integrated mixing chamber and "2 from 4" solvent selectors | 5035.0017 | Mixing Chamber Extension Kit $(\rightarrow page 136)$ |
| HPG-3400M | High-pressure gradient pump optimized for micro flows with "2 from 4" solvent selectors. The pump has no mixing chamber. | 5035.0019 | |

| Solvent Rack | Description | Part No. |
|--------------|--|-----------|
| SRD-3600 | Solvent Rack with analytical 6-channel vacuum degasser Typically for use with the following pumps: one DGP-3600 (A, AB, M, or MB) one HPG-3200 + one HPG-3400 (either A or M) in a two-stack system | 5035.9230 |
| SRD-3400 | Solvent Rack with analytical 4-channel vacuum degasser Typically for use with the following pumps: - one HPG-3400 (A or M) - two HPG-3200 (A or M) in a two-stack system. | 5035.9245 |
| SRD-3200 | Solvent Rack with analytical 2-channel vacuum degasser Typically for use with the following pumps: - one HPG-3200 (A or M - one IS0-3100A | 5035.9250 |
| SR-3000 | Solvent Rack without vacuum degasser: Typically for use with an LPG-3400 (A, AB, M, or MB) pump | 5035.9200 |

For the secure and functional positioning of the solvent reservoirs, the following Solvent Racks with integrated degasser are available:

I Tip: A Solvent Rack with an analytical degasser cannot be used with a semipreparative pump.

2.3.2 Biocompatible Pumps

The UltiMate 3000 pump series includes the following biocompatible pump versions:

- LPG-3400AB and DGP-3600AB
- LPG-3400MB and DGP-3600MB

Except for the fluidic components, the biocompatible pumps are identical to the standard pumps (stainless steel). Therefore, almost all descriptions of the standard pumps apply also to the biocompatible versions. If some detail applies to only one version, the version will be identified. The differences are as follows:

The fluidic components are made of titanium. Titanium is a base material, similar to aluminum and magnesium. When titanium is processed, a titanium oxide film builds up on the component surface, ensuring excellent corrosion resistance. Note that titanium is not as hard as stainless steel and that it has a slightly different coloration. In addition, titanium parts are lighter than parts made of stainless steel. Nevertheless, you can easily confuse titanium with stainless steel parts.

When the connection between two titanium parts is too tight, friction between the parts makes them stick together as if welded. To avoid this problem in screwed connections, Thermo Fisher Scientific uses stainless steel parts as counterparts for the titanium parts, whenever possible. Thermo Fisher Scientific recommends that you use only the capillaries shipped with the pump or original Dionex spare capillaries. The ferrules for the capillary connections are made of PEEK. Do not over tighten the connection. To ensure proper connection, the PEEK part must not deform.



Fig. 4: Installation of PEEK ferrules

Also, observe the following information:

- When substituting parts be sure to install the appropriate replacement part for the biocompatible pump version. For information about the part numbers, refer to the Service and Consumables and Spare Parts sections (→ pages 109 and 169).
- When connecting the solvent reservoirs, make sure that the filts on the end of the solvent lines are titanium frits (shipped with the pump). As standard, the accessories pack shipped with the UltiMate 3000 series Solvent Racks includes filter holders with stainless steel filter frits. Replace these frits with the titanium frits (→ page 47).
- Ring seals with a titanium spring are installed as piston seals in the biocompatible pumps, whereas common ring seals with a steel spring are used in the standard pumps. Therefore, do not confuse these seals.

2.4 Interior Components

See section 9 for the inside front panel views and fluid connections of the UltiMate 3000 pumps:

| For the | Find the | On page |
|------------------------------|--|------------|
| Isocratic pump | | |
| ISO-3100A | Interior components Fluid connections | 144 145 |
| Low-pressure gradient pumps | | |
| LPG-3400 | Interior components Fluid connections | 146 147 |
| DGP-3600 | Interior components Fluid connections | 148 149 |
| High-pressure gradient pumps | | |
| HPG-3200 | Interior components Fluid connections | 150 151 |
| HPG-3400 | Interior components Fluid connections | 153 154 |

In addition, section 9 provides information about special aspects that should be considered with a high-pressure gradient system. For information about

- Gradient combinations, see page 156
- Double Flow mode, see page 156
- the HPG-3200P semipreparative pump, see page 157

2.5 Front Panel Elements



Fig. 5: Pump front panel view

| No. | Front Panel Element | Description | |
|-----|---------------------|--|--|
| 1 | Display | Shows information about the pump, for example: General information upon power-up (→ page 59) Status screen (→ page 60) Various functions and menus (→ pages 67 and 68) Messages (→ page 96) | |
| 2 | Standby button | Switches the pump to Standby mode (the LED is red). To cancel Standby mode and resume operation, press the Standby button again (the LED is not lighted). Note : To allow the pump to change the mode, press and hold the Standby button for at least 1 second. | |
| 3 | LEDs | | |
| | Power | The LED is blue when the pump is turned on. | |
| | Connected | The LED is green when the pump is connected in Chromeleon. | |
| | Status | The LED is green when the pump is ready for operation. The LED is red when an error has been detected, for example, a leak. | |

Important: If you switch a pump to which an SRD-3x00 Solvent Rack is connected to the Standby mode, the Solvent Rack, too, will be set to Standby mode.

▲ Important: Si vous commutez une pompe à laquelle est raccordé un dégazeur SRD-3x00, en mode Veille, le dégazeur passera également en mode Veille.

2.6 Rear Panel



Fig. 6: Rear panel

| No. | Description | |
|-----|--|--|
| 1 | Power switch (\rightarrow page 19) | |
| 2 | Fuse cartridge (\rightarrow page 19) | |
| 3 | Main power receptacle (\rightarrow page 32) | |
| 4 | DC Output—Reserved for future connection of low-voltage devices. | |
| 5 | Solvent Rack port for connection of an SRD-3x00 Solvent Rack (\rightarrow page 21) | |
| 6 | Type label | |
| 7 | USB (Universal Serial Bus) port for connection to the computer for example, either direct connection or via the autosampler of the UltiMate 3000 system (\rightarrow page 19) | |
| 8 | Pressure—Analog output pressure (\rightarrow page 20) | |
| 9 | Digital I/O port for connection of an accessory, for example, a manual injection value $(\rightarrow \text{ page } 20)$ | |
| 10 | USB hub—3 additional ports for connection of one UltiMate 3000 device, such as a thermostatted column compartment of the TCC-3x00 series), or USB hub each $(\rightarrow page 19)$ | |

2.6.1 Power Switch

The power switch on the rear panel is the main power switch for the pump. Turn on the power switch before initial operation of the pump and leave it on. For routine on/off control, use the standby button on the front of the pump (\rightarrow page 17). Press and hold the button for one second to allow the pump to change the mode. Turn off the main power switch when instructed to do so, for example, before performing a service procedure or when interrupting operation for longer periods (one week or more). In this case, also observe the precautions on page 90.

2.6.2 Fuse Cartridge

The fuse cartridge contains two slow-blow fuses rated at 2 A, 250 V. For information about how to change the fuses, see page 131.

2.6.3 USB Connector

The Chromeleon Chromatography Management System can use a USB connection to control the pump. Data is transferred digitally via the appropriate USB cable (\rightarrow page 32). To ensure trouble-free operation, use only the cables shipped with the pump.

The internal USB hub (\rightarrow Fig. 6, no. 10) allows you to connect three other instruments in the UltiMate 3000 product line or three external USB hubs to the pump.

- ▲ Important: Thermo Fisher Scientific recommends using these USB ports for connections to Dionex instruments only. Thermo Fisher Scientific cannot guarantee correct functioning if instruments from other manufacturers are connected.
- Important: Thermo Fisher Scientific recommande d'utiliser les ports USB uniquement pour les raccordements aux instruments Dionex. Thermo Fisher Scientific ne peut garantir le bon fonctionnement si les instruments d'autres fabricants sont raccordés.

For information about how to connect the pump to the Chromeleon computer, see sections 3.4.1 and 3.4.2 (\rightarrow page 32).

2.6.4 Pressure (Analog Pressure Output)

The analog pressure output records the operating pressure of the pump. The pressure output is set to 5 mV/bar (50 mV/Mpa). You may connect a device, such as the UCI-100 Universal Chromatography Interface, a recorder, or an A/D converter, to monitor the pump pressure.

If a DGP-3600 is operated by Chromeleon, use the **AnalogOut** property to determine whether the pressure of the right or left pump is available at the analog pressure output. (Chromeleon does not support this property for the other pumps in the UltiMate 3000 pump series.)

| Pin Assignment for 2-Pin Cinch Connector | | |
|--|-------------------|--|
| Inner ring: | Signal (pressure) | |
| Outer ring: | GND | |

The analog pressure output always records the pressure of the pump. This is also true when the pump is operated together with am FLM-3x00 Flow Manager in an UltiMate 3000 system. This means that the pump pressure is recorded (not the column pressure).

2.6.5 Digital I/O

The digital I/O port provides three inputs and four relay outputs that can be used to trigger or read in external events.

You can use the inputs (1-3) as universal inputs and read them in Chromeleon. The outputs (relays 1-4) can be used as universal outputs, they can be controlled via Chromeleon, or they can be assigned pump-internal special functions. For more information, see section 5.6.4 (\rightarrow page 88).

Important: The maximum switching voltage of the relays is 24 V. The switching current must not exceed 100 mA.

Important: La tension maximale de commutation des relais est de 24 V. L'intensité de commutation ne doit pas dépasser 100 mA.

For information about the pin assignment and the signal levels, see page 186.

2.6.6 Solvent Rack

Use this port to connect an SRD-3x00 Solvent Rack with integrated analytical vacuum degasser to the pump. A Solvent Rack with an analytical degasser cannot be used with the HPG-3200P semipreparative pump

- **Important:** Do not substitute any other Solvent Rack for the Solvent Racks mentioned in the table on page 14.
- **Important:** Ne remplacez les dégazeurs de la série SRD-3x00 mentionnés sur la page 14 par aucun autre type de dégazeur.

For information about the pin assignment of the Solvent Rack port, see page 183. For information about how to install and operate the Solvent Rack, see the *Operating Instructions* for the instrument.

2.7 Fluid Connections

The front panel tilts upward to provide easy access to the fluid connections in the pump. The open cover locks in the topmost position. For an overview of the inside front panel views and the fluid connections of the UltiMate 3000 system series pumps, see page 143 and the following pages.

Important: Do not place any heavy objects on the open front panel door. This may damage the door.
 When lifting or moving the pump, always lift by the bottom or sides of the instrument. Lifting the pump by the front panel may damage the front panel door.
 Important: Ne placez aucun objet lourd sur la porte ouverte du panneau avant. Ceci peut endommager la porte.
 Lorsque vous soulevez ou déplacez la pompe, saisissez la toujours par le dessous ou les côtés de l'instrument. Soulever la pompe par le panneau avant risque d'endommager la porte du panneau avant.

For information about the connections in the low-pressure and high-pressure sections of the pump, see page 183.

2.8 Rear Seal Wash System

The pump is equipped with an active real seal wash system. When using buffer solutions, Thermo Fisher Scientific recommends continuously rinsing the back of the piston seal to remove salt crystals and prolong the life of the seal.

For the standard setup, the real seal wash system consists of a peristaltic pump and an internal liquid reservoir with integrated detector. Fig. 7 shows how the components are connected upon shipment of the pump.



Fig. 7: Rear seal wash system (standard setup) Top: ISO-3100A and LPG-3400; bottom: all other pumps types (here DGP-3600)

As an alternative to the standard setup, rear seal washing can be performed also from an external liquid reservoir. For installation instructions, see page 52.



Fig. 8: Rear seal wash system (alternative setup), here for a DGP-3600

Use this setup, for example, if you

- want to use fresh wash liquid for each wash cycle.
- use solvents with a high salt content.
- observe desposits on the pistons or increased piston seal wear with specific solvents.

2.9 Outlet Unit

The outlet unit comprises a pressure transducer for the system pressure, purge valve, and inline filter (filter holder with filter frit). It depends on the pump type whether the outlet unit includes a mixing chamber (\rightarrow page 24).

In a DGP-3600, each of the two pumps is fitted with a separate outlet unit. Do not interconnect the outlet units. Always direct them to separate fluid systems.



Fig. 9: Outlet unit (here with mixing chamber)

2.9.1 Inline Filter

The pump is shipped with the appropriate filter frits, depending on the pump with a porosity of 0.5 μm or 10 μm

LPG-3400M, LPG-3400MB, DGP-3600M, and DGP-3600 MB

These pumps are mainly used in nano and capillary HPLC applications, together with an FLM-3x00 Flow Manager. To protect the thin capillaries, the flow splitter, and flow control valve in the flow manager, these pumps have filter frits with a porosity of $0.5 \,\mu$ m installed.

All other pumps

All other pumps have filter frits with a porosity of 10 μ m installed.

| Filter Frits with a Porosity of | Part No. |
|--|-----------|
| 0,5 µm for LPG-3400M and DGP-3600M | 6000.0045 |
| 2 µm for LPG-3400MB and DGP-3600MB | 6268.0036 |
| 10 μ m for LPG-3400AB and DGP-3600AB | 6268.0032 |
| 10 μm for all other pumps | 6268.0031 |

The accessories kit for some pump types includes filter frits with different porosities. Therefore, make sure not to confuse the different frit types.

2.9.2 Mixing Chamber

The following pumps are equipped with a dynamic mixing chamber (volume: 328 μ L) in the outlet unit (\rightarrow Fig. 59, page 135)

- LPG-3400A and LPG-3400AB
- DGP-3600A and DGP-3600AB (These pumps have two mixing chambers.)
- HPG-3x00A

For poorly miscible solvents, higher flow rates, or special applications, you can easily extend the mixing chamber volume (by 600 μ L or 1200 μ L) by installing the associated extension kit that is available as an option (\rightarrow pages 135 and 136).

LPG-3400A and DGP-3600A only

For gradient separations at low flow rates (for example, less than 100 μ L/min) or very steep gradients, you can operate these pumps without mixing chamber (\rightarrow page 138), and thus adapt the gradient delay volume to your requirements. Reducing the gradient delay volume allows changes in the solvent composition to become effective on the column much earlier, thus having a faster effect on the elution of analytes.

The gradient delay volume of an HPLC system is defined as the volume that the pump must deliver until a change in solvent composition reaches the head of the column. For information about the delay volume of the different pumps, see page 158 and the following pages.

HPG-3200P only

As a standard, this pump is shipped with an 328 μ L mixing chamber plus 600 μ L extension. An 1200 μ L mixing chamber extension is available as an option (\rightarrow page 136).

HPG-3x00M only

A static mixer is available from Thermo Fisher Scientific as an option for the HPG-3x00M pumps (\rightarrow page 141). The mixer improves the mixing quality of the 2 combined solvent streams delivered by the pump, and thus guarantees a smoother baseline.

2.10 Leak Sensor

A leak sensor is installed inside the pump. If liquid collects in the drip tray under the fluid connections, the leak sensor reports a leak, and the **Status** LED on the front panel door changes to red. It depends on the leak sensor mode setting (\rightarrow page 84) whether a message appears on the front panel display and a beep alerts you.

When the leak sensor reports a leak, eliminate the cause for the leakage and dry the leak sensor (\rightarrow page 110). To remove the message from the display, select **Clear** on the navigation bar (\rightarrow page 70).

2.11 Vacuum Degasser

Usually, a vacuum degasser is used to remove gas bubbles trapped in the solvents. With the LPG-3400 pump models, the vacuum degasser is integrated in the pump. For the DGP-3600 pump models, the HPG-3x00A and M pumps, and the ISO-3100A pump, Thermo Fisher Scientific recommends using the appropriate SRD-3x00 Solvent Rack with integrated degasser (\rightarrow page 14) or any other external vacuum degasser.

Normal phase eluents usually show only a low concentration of dissolved gases. Therefore, it is normally not required to use a degasser with these eluents.

I Tip: The degassers run quietly. Even if the vacuum pump is running, the operating noise is very low. Higher frequencies may slightly increase the noise level, but this does not impair the degassing performance.

When using a HPG-3200P pump, make sure to connect a preparative degasser (not an analytical degasser).

2.12 Chromeleon Software

The pump can be operated with the Chromeleon Chromatography Management System. To control the pump, an appropriate Chromeleon version and a **Timebase Class 1** Chromeleon license are required. Two modes of software control are available:

• Direct Control

With direct control, you select operating parameters and commands in the **Commands** (F8) dialog box or on a control panel. Direct commands are executed as soon as they are entered. For routine operation, most parameters and commands are available also on a control panel. For more information about direct control, see page 62.

Automated Control

With automated control, you create a program (or PGM File). This is a list of control commands, executed in chronological order, for automated operation of the pump. Programs can be created automatically with the help of a software wizard or manually by editing an existing program. For more information about automated control, see page 65.

I Tip: All software details in this manual refer to *Chromeleon 6.80*.

If you want to operate the pump with Chromeleon 7, refer to the following documents for information about how to perform the related processes in Chromeleon 7 (all documents are included in the Chromeleon 7 shipment):

- *Chromeleon 7 Help*—provides extensive information and comprehensive reference material for all aspects of the software.
- *Quick Start Guide*—describes the main elements of the user interface and guides you step-by-step through the most important workflows.
- *Reference Card*—provides a concise overview of the most important workflows.
- *Installation Guide*—provides basic information about module installation and configuration. For specific information about a certain module, refer to the *Chromeleon 7 Instrument Configuration Manager Help*.

Note the following:

- Chromeleon 7 terminology is different from the terminology used in Chromeleon 6.80. For details, refer to the 'Glossary Chromeleon 7,' which is available in the Documents folder of your Chromeleon 7 installation.
- Chromeleon 7 may not yet support all functions supported in Chromeleon 6.80.
2.13 System Wellness, Predictive Performance, and Diagnostics

System Wellness monitors the health of the pump. Therefore, the pump supports several performance and reliability features that can help you detect small problems before they turn into big ones:

- Internal monitoring of all mechanical operations
- Automatic self test upon power up
- Leak sensor (\rightarrow page 25)
- Active rear seal wash system (\rightarrow page 22)
- Monitoring of the liquid level for rear seal washing (\rightarrow Case B, page 49)

When an error is detected, the **Status** LED on the front panel changes to red and a message appears on the front panel display (\rightarrow page 96).

When the pump is operated with Chromeleon, additional functions for estimating the lifetime of consumables and monitoring and recording service and (re)qualification information (= predictive performance) are available (\rightarrow page 86). To check the performance of certain pump components and the overall performance of the instrument, Chromeleon also supports diagnostic functions for the pumps of the UltiMate 3000 system (\rightarrow page 87).

3 Installation

3.1 Facility Requirements

- Make sure that the installation site meets the power and environmental specifications listed in the Technical Information section (→ page 159).
- Install the pump in the laboratory on a stable surface that is free of vibrations.
- Make sure that the surface is resistant to solvents.
- Avoid locations with extreme changes in temperature (such as direct sunlight or drafts) and high humidity.
- Allow sufficient clearance behind and to the sides of the pump for power connections and ventilation.

3.2 Unpacking

All electrical and mechanical components of the pump are carefully tested before the instrument is shipped from the factory. After unpacking, please inspect the instrument for any signs of mechanical damage, which might have occurred during transit.

1 Tips: Immediately report any shipping damage to both, the incoming carrier and Thermo Fisher Scientific. Shipping insurance will compensate for the damage only if reported immediately.

Keep the original shipping container and packing material. They provide excellent protection for the instrument in case of future transit. Shipping the unit in any other packaging automatically voids the product warranty.

- 1. Place the shipping container on the floor and remove the accessories kit and the power cord.
- 2. Grasp the pump by the sides. Slowly and carefully, pull the instrument out of the shipping container and place it on a stable surface.
 - **M** Important: To prevent the pump from falling, always lift by the bottom or sides of the unit. Do not lift the unit by the packaging material or the front panel door.
 - ▲ Important: Afin d'empêcher la pompe de tomber, saisissez-la par le bas ou les côtés. Ne soulevez la pompe à l'aide du matériau d'emballage ou par la porte du panneau avant.

- 3. Remove the foam spacers, and then remove the polythene packaging.
- 4. Tilt the front panel upward and remove the foam inserts securing the front panel door during shipment.
- 5. Before connecting the pump to the power source, wait approximately 4 hours to allow the instrument to come to room temperature and to allow any condensation that might have occurred during shipping to evaporate. After 4 hours, check the pump; if condensation still exists, allow the pump to continue to warm up (without connecting it to the power source) until the condensation is completely gone.

3.3 Positioning the Pump in the UltiMate 3000 System

If the pump is part of an UltiMate 3000 system, for example, for analytical HPLC applications, Thermo Fisher Scientific recommends stacking the individual modules, for example, as shown in Fig. 10 and interconnecting them on the rear panel as shown in Fig. 11.

However, the arrangement of the system modules depends on the application. If the pump is part of an UltiMate 3000 Proteomics MDLC system for nano or capillary HPLC, the UltiMate 3000 Proteomics MDLC system manual provides information about how to arrange and connect the modules for these applications.



Fig. 10: Example for an UltiMate 3000 system



Fig. 11: Example for the rear panel connections on an UltiMate 3000 system (here with VWD-3x00)

1 Tip: Apart from the Solvent Rack, all modules of the UltiMate 3000 system can be connected separately to the computer. However, Thermo Fisher Scientific recommends interconnecting all modules, and then connecting the system to the Chromeleon computer via only one connection. To do so, use the USB hub on the pump or autosampler.

3.4 Connecting the Pump

3.4.1 General Information

If you want to operate the pump with Chromeleon

Verify that Chromeleon is installed on the computer and that the license code is entered *before* you connect the pump to the USB port on the Chromeleon computer and turn on the pump power. Only if you install Chromeleon first, the USB driver for the pump is automatically loaded and the Windows[®] operating system can detect the pump when the power is turned on.

3.4.2 Connecting the USB Cable

Connect the pump to the Chromeleon computer via the USB ports on the rear panel (\rightarrow Fig. 6, page 18). To do so, select one of the following alternatives:

- Connect the pump directly to the USB port on the computer.
- Connect the pump to an internal USB port on another module in the UltiMate 3000 system that is connected to the computer.
- Connect the pump to the computer via an external USB hub.

The following cables are available:

| Cable | Part No. |
|-----------------------------------|-----------------------------------|
| USB cable, 5 m, A to type B cable | 6911.0002 (shipped with the pump) |
| USB cable, 1 m, A to type B cable | 6035.9035 |

Tip: The USB standard limits the USB cable length to 5 meters. Each USB device can be separated from the PC or next USB hub by no more than 5 meters.

3.4.3 Connecting the Power Cord

Use the power cord shipped with the pump to connect the instrument to the main power source. Connect the power cord from the main power receptacle on the rear panel (\rightarrow Fig. 6, page 18) to the power source that is connected to a true ground. No manual adjustment is required to adapt the line voltage to local voltage requirements.

3.4.4 Connecting the Solvent Rack and Digital I/O

Solvent Rack

If the UltiMate 3000 system includes an SRD-3x00 Solvent Rack, connect this port on the rear panel of the pump to the related port on the rear panel of the Solvent Rack. The appropriate connection cable is included in the accessories pack of the Solvent Rack.

Digital I/O

If you want to use the inputs and relay outputs in Chromeleon (\rightarrow page 88), connect the Digital I/O port on the rear panel of the pump to the associated port on the Chromeleon computer.

3.5 Setting Up the Pump in Chromeleon

This section provides brief instructions for setting up Chromeleon. For details about any of these steps, see the Chromeleon Help.

Tip: When the pump is connected to the Chromeleon computer, verify that the Chromeleon software is installed *before* turning on the pump power for the first time. Only then, the Windows operating system will detect the pump automatically when the pump power is turned on.

3.5.1 Loading the USB Driver for the Pump

- 1. Turn on the computer power, if it is not already on.
- 2. Under Windows Vista[®] (Windows[®] XP, Windows[®] 7, or Windows[®] Server 2008) log on as a
 - Local administrator if the computer is a local computer.
 - User with local computer administrator privileges if the computer is a network computer.
- Start the Chromeleon Server Monitor program by double-clicking the Chromeleon Server Monitor icon ion on the Windows taskbar.

If the Server Monitor icon is not on the taskbar, click **Start** on the taskbar, point to **Programs** (or **All Programs**, depending on the operating system), point to **Chromeleon**, and then click **Server Monitor**.

- 4. Click **Start** to start the server.
- 5. Click **Close** to close the Server Monitor window. The Server Monitor icon appears on the taskbar.



Tip: Clicking the **Quit Monitor** button quits (exits) the Server Monitor program, but does not stop the server. To stop the server, click **Stop**.

6. Turn on the main power switch on the rear panel of the autosampler.

7. *Windows Vista, Windows 7, and Windows Server 2008* will automatically detect the new pump and perform the USB installation.

If Windows Vista fails to detect the pump and launches a wizard instead, this indicates that you connected the pump to the computer and turned on the power for the first time *before* you installed Chromeleon. To resolve the problem:

- a) Click **Cancel** to exit the wizard.
- b) Turn off the pump and unplug the USB cable from the computer.

- c) Install Chromeleon.
- d) Reconnect the USB cable to the computer and turn on the power to the pump. Vista will now detect the pump and install the USB software for the pump automatically.
- 8. Windows XP

will automatically detect the new pump and launch the **Found New Hardware Wizard**, which guides you through the USB installation. Select the following options:

- a) If asked whether Windows can connect to Windows Update to search for software, select **No, not this time**.
- b) Accept the default option (Install the software automatically) and click Next>.
- c) Click **Finish** when the wizard reports that the software for the pump has been installed.

If Windows fails to detect the pump and a message box asks for a USB configuration file (cmwdmusb.inf), this indicates that you connected the pump to the computer and turned on the power for the first time *before* you installed Chromeleon.

To resolve the problem:

- a) Click Cancel in the Windows message box.
- b) Turn off the pump and unplug the USB cable from the computer.
- c) Install Chromeleon.
- d) Reconnect the USB cable to the computer and turn on the power to the pump. Windows will now automatically detect the pump and launch the **Found New Hardware Wizard.**

3.5.2 Installing the Pump

After the USB software for the Pump has been installed (\rightarrow page 34), install and configure the pump in Chromeleon:

- 1. Start the Chromeleon Server Monitor if it is not already running (\rightarrow page 34)
- 2. Start the Chromeleon Server Configuration program by clicking Start on the taskbar. Point to Programs (or All Programs, depending on the operating system), point to Chromeleon, and then click Server Configuration.
- 3. If necessary, click the plus sign next to the server name to display the items underneath.
- 4. Select the timebase to which the pump will be assigned, or create a new timebase (on the **Edit** menu, click **Add Timebase**).
- 5. Open the **Add device to timebase** dialog box. To do so, click **Add Device** on the **Edit** menu or right-click the timebase and click **Add Device** on the menu.
- 6. On the **Manufacturers** list, click **Dionex HPLC: UltiMate 3000** and on the **Devices** list, click the pump: **DGP-3600 Pump**, **HPG-3x00 Pump**, **ISO-3100 Pump**, or **LPG-3x00 Pump**.
- 7. A wizard guides you through the installation. Chromeleon connects to the pump and transfers the settings from the instrument firmware to Chromeleon, setting the options on the wizard pages accordingly.
- 8. On each wizard page, verify that the settings are correct and select additional settings if needed. For a description of the wizard pages, see section $3.5.3.1 (\rightarrow \text{ page } 37)$.
- 9. Click **Finish** to complete the installation of the pump.
- 10. On the **File** menu, click **Save Installation** and then close the Server Configuration program.

3.5.3 Configuring the Pump

3.5.3.1 Initial Installation

During the installation, Chromeleon connects to the pump and transfers the settings from the instrument firmware to Chromeleon, setting the options on the wizard pages accordingly. Verify that the default settings are correct and make additional settings if needed. You may reopen the configuration pages later again to change the settings (\rightarrow page 44).

Tip: Changing the settings for a specific application in the **Commands** dialog box, in a program file (PGM), or on a control panel will not change the default settings on the configuration pages.

For additional information about a page, click Help.

General Page

Define the general instrument parameters.

| General | | × |
|-------------------------|--------------------------------------|-----------------|
| | | |
| | | |
| | | |
| | | |
| | Firmware Download | |
| | | |
| Module <u>A</u> ddress: | | |
| | | B <u>r</u> owse |
| | | |
| | | |
| | | |
| | | |
| | | |
| | < <u>B</u> ack <u>N</u> ext > Cancel | Help |
| | | |

Fig. 12: General page

• Demo Mode

Verify that the check box is cleared. In the Demo Mode, Chromeleon simulates the functions of the pump. If the Demo Mode is enabled, the **Module Address** input field will be unavailable. If you exit this page without having entered a module address, the Demo Mode will be enabled automatically.

Module Address

To enter the module address of the pump, click **Browse** and then double-click the pump that you want to use. The address is automatically entered in the **Module Address** field.

• Firmware Download

Clicking this button updates the pump firmware with the version available in Chromeleon. (The button appears dimmed if the Demo Mode is enabled.)

A message displays the version of the currently installed firmware version and of the version that will be downloaded from Chromeleon. Click **OK** to start the download.

For the download, make sure that

- the Chromeleon server is in **running idle** mode.
- the pump is unpressurized (for example, the purge valve on the outlet unit is open) and that the flow is turned off.

The download, which can take several minutes, is complete when **Download finished successfully** appears in the **Messages Server** window in the Chromeleon Server Configuration program. (The message appears also in the Chromeleon Audit Trail.)

Immediately after the new firmware has been downloaded from Chromeleon to the instrument, the pump performs a reset. For about 15 seconds, the internal boot loader is updated. Therefore, do not turn off the pump while the boot loader is updating

| ▲ Important: | During the download, make sure that the communication between the pump and Chromeleon is not interrupted and do not turn off the instrument. This may lead to malfunction of the pump. If the download is not successful, the related messages appear in the Audit Trail. In this case, turn off the pump. Turn on the pump again and repeat the download as described above. If the download fails, contact Thermo Fisher Scientific Service for Dionex HPLC Products. |
|--------------|--|
| ▲ Important: | Au cours du téléchargement, assurez-vous que la communication entre la pompe et Chromeleon n'est pas interrompue et n'arrêtez pas l'instrument. Ceci peut entraîner des dysfonctionnements de l'instrument. Si le téléchargement échoue, arrêtez l'instrument. Remettez l'instrument en marche et répétez l'opération de téléchargement tel que décrit ci-dessus. Si le téléchargement échoue, contactez le service après-vente Thermo Fisher Scientific Service pour Dionex HPLC Produits. |

When you leave the **General** page during initial installation of the pump (with the Demo Mode being disabled), Chromeleon attempts to connect to the pump, retrieve the settings from the pump firmware, and set the options (pump type and pump head) on the corresponding wizard pages. Confirm the related message with **OK**.

Devices Page

| evices | | |
|---------------------|-----------------------------|------------------------------|
| | | |
| Device <u>N</u> ame | e: Pump | |
| Left Pump: | LoadingPump | on ComputerName_UltiMate_1 |
| | uses FLM-3x00 Flow Splitter | <none></none> |
| | purges via WPS-3000SL | UM3WPS_PURGE0 |
| Right Pump: | MicroPump | on ComputerName_UltiMate_2 💌 |
| | uses FLM-3x00 Flow Splitter | FLM3x00_FLOWSPLITTER_1 |
| | purges via WPS-3000SL | <none></none> |
| 🔽 Share <u>e</u> lu | ient bottles | |
| 🔽 Share <u>w</u> a | aste bottle | UM3WPS_PURGE1 |
| ✓ Pressure | Signal(s) | |
| | | |
| | | |
| | < <u>B</u> ack <u>N</u> ext | t> Cancel Help |

The pump type determines which settings can be made on this page.

Fig. 13: Devices page (here: DGP-3600M)

- All pumps
 - Device Name

Displays the name under which the pump is identified in the installation environment and in the Chromeleon client. The default name is **Pump** (**MicroPump** for an LPG-3400M(B)). To control the pump with the existing control panels, accept the default name. If you enter a different name, you may have to re-link the controls on the control panels and edit the device name in the program files.

purges via WPS-3000SL¹

If the UltiMate 3000 system includes a WPS-3000SL autosampler, use the field to specify whether the pump is purged via the membrane pump of the autosampler. The default setting is **<None>**.

All pumps except DGP-3600

Select UM3WPS_PURGE0 to connect the pump fluidically with the autosampler.

¹ For patent reasons, this function must not be used in the following countries: China, Germany, Great Britain, Japan, and the USA.

DGP-3600 only

To fluidically connect the two pumps of a DGP-3600 alternately with the autosampler, select UM3WPS_PURGE0 for one pump and UM3WPS_PURGE1 for the other pump. To connect only one pump, select UM3WPS_PURGE0 for the pump to connect and select <None> for the other pump.

For more information about how to purge the pump via the autosampler, see page 56.

Pressure Signal(s)

The check box is selected by default. Accept this setting if you want to record the pump pressure. Chromeleon generates the appropriate channels for data acquisition. For more information, see section 5.5.6 (\rightarrow page 80).

• DGP-3600 only

• Left Pump and Right Pump

Display the names under which the two pumps of a DGP-3600 are identified in the installation environment and in the Chromeleon client.

For the DGP-3600M(B), the default names are **LoadingPump** for the left pump and **MicroPump** for the right pump. For the DGP-3600A, the default names are **PumpLeft** and **PumpRight**. To control the pumps with the existing control panels, accept the default names. If you enter different names, you may have to re-link the controls on the control panels and edit the device names in the program files.

♦ on

Select the timebase to which the related the pump is assigned. The default setting is the timebase in which the pump is installed. To change the setting, click the arrow next to the on field and select a different timebase from the drop-down list. The list shows all timebases for that server.

Share eluent bottles

The check box is selected by default. Accept this setting if both pumps of the DGP-3600 are connected to the same set of solvent reservoirs. Clear the check box if the pumps are connected to different sets of reservoirs. In this case, Chromeleon supports the %A/B/C_RemainTime and %A/B/C_WarningLimit properties separately for each pump. If the check box is selected, Chromeleon assigns the properties the same values for each pump.

• Share waste bottles

The check box is selected by default. Accept this setting if both pumps are connected to the same waste container. Clear the check box if the pumps are connected to different waste containers. In this case, Chromeleon supports the

Waste_RemainTime and Waste_WarningLimit properties separately for each pump. If the check box isselected, Chromeleon assigns the properties the same values for each pump

• DGP-3600M(B), LPG-3400M(B), HPG-3200(A and M) and HPG-3400 only

uses FLM-3x00 Flow Splitter

If the UltiMate 3000 system includes an FLM-3x00 Flow Manager, use these fields to specify whether the pump is connected to a flow splitter. The default entry is **<None>**. To change the setting, click the arrow next to the input field and select the splitter from the drop-down list.

When you assign a flow splitter to a pump, the pump flow settings apply directly to the flow on the splitter outlet, that is, the split ratio is considered automatically. In the standard configuration of the UltiMate 3000 system with a DGP-3600M(B) pump, Thermo Fisher Scientific recommends assigning the flow splitter to the right pump (MicroPump).

Configuration complete Page

Basic configuration of the pump is complete. You can finish the installation of the pump and or continue with the wizard for the more advanced settings:

- The default setting is Accept all settings and close the configuration wizard. Click Finish if you want to complete the configuration and accept the default values for the more advanced settings. Afterward, click Save on the File menu and then close the Server Configuration program.
- Select **Continue with wizard to configure advanced settings** to continue with the advanced settings (flow limits, pressure limits, and solvent names) and determine which relays and digital inputs shall be available in Chromeleon. See the following sections for a description of these settings.

Head Type & Limits Page or Right Pump and Left Pump Pages

Check and change if necessary the upper and lower limits for the flow rate, the pressure unit, and the pressure limits.

The pump type determines which pages are available. These are **Right Pump** and **Left Pump** for a DGP-3600 and **Head Type & Limits** for all other pump types.

| Minimum: 0.000 ml/min Maximum: 10.000 ml/min essure Image 0500 bar Unit: bar Minimum: 0 bar Image 0500 bar Maximum: 400 bar Image 0500 bar | 1010 | Range 0.000 |) 10.000 ml/min | | uble flow | |
|--|-------|-------------------|-----------------|---------------|-----------|--|
| Maximum: 10.000 ml/min essure Range 0 500 bar Unit: bar Minimum: 0 bar Maximum: 400 bar | ļ | <u>M</u> inimum: | 0.000 | ml/min | | |
| essure Range 0 500 bar Unit: bar 💌 Minimum: 0 bar Maximum: 400 bar | | M <u>a</u> ximum: | 10.000 | ml/min | | |
| Range 0 500 bar Unit: bar 💌 Minjimum: 0 bar Maximum: 400 bar | ressu | re | | | | |
| Mi <u>n</u> imum: 0 bar Ma <u>x</u> imum: 400 bar | | Range () 51 | 00 bar | <u>U</u> nit: | bar | |
| Ma <u>x</u> imum: 400 bar | | Mi <u>n</u> imum: | 0 | bar | | |
| | | Ma <u>x</u> imum: | 400 | bar | | |

Fig. 14: Head Type & Limits page (here: HPG-3x00A)

• All pumps

The box on the top left indicates the pump version (installed head type): Analytical, Micro, or Semipreparative. The head type affects the maximum allowed flow rate and pressure.

♦ Flow

Shows the allowed flow rate range. Under **Minimum** or **Maximum**, enter a new value to change the lower or upper limit for the flow rate range.

◆ Pressure

Shows the allowed pressure range. Under **Unit**, change the pressure unit (psi, bar or MPa) if required. Under **Minimum** or **Maximum**, enter a new value to change the lower or upper limit for the pressure range.

• HPG-3x00 only

The **High pressure gradient** check box is selected by default. Only then, **Double flow** mode can be enabled. In Double flow mode, both pump blocks are used for the delivery. In this way, the maximum allowed flow rate is doubled (\rightarrow page 156).

Solvents Page or Right Solvents and Left Solvents Pages

Check and change if necessary the number of solvents delivered by the pump and the solvent names. The pump type determines which pages are available. These are **Right Solvents** and **Left Solvents** for a DGP-3600 and **Solvents** for all other pump types.

• Number of Solvents

Indicate how many solvents are delivered by the pump. The pump type determines how many solvents can be delivered.

| Pump | Maximum Number of Solvents |
|-----------------------|-------------------------------|
| ISO-3100 | 1 |
| HPG-3200 | 2 |
| LPG-3400 and HPG-3400 | 4 |
| DGP-3600 | 2x3 |

• Solvent Names

Enter a name for each connected solvent. A maximum of 30 characters is permitted. The names of the solvents appear, for example, in the gradient display of the online control panel and in the report.

Relays Page

This page lists all available relays. Select a check box to enable the corresponding relay. If a check box is cleared, the relay will not be available in Chromeleon. To change the relay name or timebase assignment, double-click the relay name (or press the F2 button) and make the changes in the **Devices Configuration** dialog box. For more information about the relays, see page 20.

Inputs Page

This page lists all available remote inputs. Select a check box to enable or disable the corresponding remote input. If a check box is cleared, the input will not be available in Chromeleon. To change the input name or timebase assignment, double-click the input name (or press the F2 button) and make the changes in the **Devices Configuration** dialog box.

3.5.3.2 Changing the Configuration Properties

To change the default configuration settings, reopen the configuration pages:

- 1. Start the **Server Configuration** program (\rightarrow page 36).
- 2. Right-click the pump in the timebase and select **Properties** on the menu.
- 3. Change the settings as needed. For a description of the wizard pages, see section 3.5.3.1 (→ page 37).

On the **General** page, a **Retrieve configuration from module** button is available. Click this button to transfer the current pump configuration to Chromeleon. This is necessary only if you connect a different pump or if automatic matching was not performed correctly when the wizard was started.

In addition to the pages provided during the initial installation of the autosampler, Chromeleon provides an **Error Levels** page. This page classifies the severity of any errors that occur. It is generally not necessary to change the default settings.

4. To save the changed configuration, click **Save** on the **File** menu and then close the **Server Configuration** program.

4 Preparation for Operation (Startup)

4.1 Overview of Actions

- Important: The pump is filled with 2-propanol when being shipped from the factory. During initial operation of the pump, make sure that the solvents used are miscible. Otherwise, use an appropriate intermediate solvent.
- Important: La pompe est stockée sous 2-propanol lorsqu'elle est expédiée depuis l'usine. Lors du démarrage initial de la pompe, assurez-vous que les solvants utilisés sont miscibles. Dans le cas contraire, utilisez un solvant intermédiaire approprié. Même d'infimes particules peuvent endommager le système.

After you have unpacked, positioned and connected the pump as described in sections 3.1 through 3.4 (\rightarrow page 29 and following pages), prepare the pump for operation. Follow the sequence of steps below:

1. The fluid connections depend on your application. Connect the pump to the autosampler and to the column compartment or flow manager. If the UltiMate 3000 system does not include an autosampler, you may as well connect a manual injection valve to the pump (\rightarrow page 184).

When connecting capillaries to the pump, observe the general precautions on page 46

If you want to connect the pump to the flow splitter of the flow manager, use the special capillary from the respective application kit. Do not substitute this capillary for any other capillary. For more information, see the *Flow Manager Manual*.

- 2. Install the Solvent Rack if applicable. For installation details, see the *Solvent Rack Manual*.
- 3. Connect the solvent reservoirs, and then connect the solvent lines to the degassing channels (\rightarrow page 47).
- 4. Verify that the peristaltic tubing is engaged in the peristaltic pump and fill the liquid reservoir of the rear seal wash system (\rightarrow page 49).
- 5. Connect drain tubing for discharging liquid leaks that might have accumulated in the interior to the drain port at the bottom right of the pump (\rightarrow page 53).
- 6. If you want to operate the pump with Chromeleon Set up the pump in Chromeleon if it is not already set up (\rightarrow page 34).
- 7. Turn on the pump (\rightarrow page 60).
- 8. Turn on rear seal washing (\rightarrow page 79).

- 9. Purge the pump (\rightarrow page 54).
- 10. Check and change the leak sensor setting if necessary (\rightarrow page 84).
- 11. Adjust the brightness and contrast of the front panel display if necessary (\rightarrow page 85).
- 12. Before using the pump for sample analysis, equilibrate the entire system (\rightarrow page 57).

4.2 General Precautions for Connecting Capillaries

When connecting capillaries to the pump, observe the following general precautions:

- When connecting capillaries, make sure that the connectors are free from contaminants. Even minute particles may cause damage to the system
- Capillary outlets, which may be capped by black rubber stoppers, are provided in the enclosure bottom. Through these rubber stoppers and outlets, direct the capillary from the outlet unit of the pump to the outside.



Fig. 15: Capillary outlets

Slit the stopper if necessary and thread the capillary through the slot.

- Reuse used fittings and ferrules only for the same capillary connection. This is to avoid increased dead volume.
- Different fitting systems are used in an UltiMate 3000 system. Therefore, install the capillaries and fittings only at the positions for which they are intended.
- Use only the capillaries shipped with the pump and original Dionex spare capillaries.

4.3 Solvent Reservoirs

For the secure and functional positioning of the solvent reservoirs, the UltiMate 3000 system series includes Solvent Racks with and without integrated vacuum degasser (\rightarrow page 14). All Solvent Racks are shipped with solvent reservoirs and appropriate tubing, including frit holders with filter frits.



Fig. 16: Pump with Solvent Rack

If the UltiMate 3000 system includes an UltiMate 3000 series autosampler, Thermo Fisher Scientific recommends degassing the wash liquid on a continuous basis, via the degasser of either the LPG-3400 pump or an appropriate SRD-3x00 Solvent Rack. The procedure how to prepare and install the wash liquid lines is similar to the steps below. For more information, see the *Autosampler Manual*.

4.3.1 General Precautions

When connecting the solvent reservoirs, observe the following general precautions:

- Always install filter frits on the solvent lines. This prevents contaminants from reaching the HPLC system.
- Regularly check the filter frits for permeability. Replace the filter frits at regular intervals. This is especially important when using aqueous solvents. Aqueous solvents may contaminate the filters with algae and other microorganisms that deposit on the filter frits. Therefore, also replace the solvents at regular intervals. Rinse the reservoirs thoroughly before refilling them. Replace the filter frits as necessary.
- As standard, the filter frits shipped with the Solvent Rack are 10 µm stainless steel frits. If the UltiMate 3000 system includes a biocompatible pump (→ page 14), replace the stainless frits with the titanium frits from the accessories kit of the pump. To do so, unscrew the top part of the filter holder from its bottom part and replace the filter frit. When placing the new filter frit into the bottom part, make sure that the frit is in a level position (avoid tilting the frit).

The same applies if the system includes a semipreparative HPG-3200P pump. In this case, too, use the special frits from the accessories kit.

- Make sure that the tubing connecting the pump to the degasser is as short as possible and locate the solvent reservoirs as close as possible to the pump. To avoid formation of gas bubbles in the reservoirs and reformation of gas bubbles in the solvent, make sure that the reservoirs are on the same level or higher as the pump. Therefore, stack the rack onto the pump as shown in Fig. 16 (→ page 47).
- Normal phase eluents usually show only a low concentration of dissolved gases. Therefore, it is normally not required to use a degasser with these eluents.

4.3.2 Connecting the Solvent Reservoirs

- 1. Feed the solvent line through the retaining guide, which holds the tubing in place in the bottle, and then into the open hole in the reservoir cap.
- 2. Slide the filter holder with filter frit onto the end of the solvent line. Verify that the end of the solvent line is cut straight and is not deformed. Cut the tubing straight if necessary. Use only the original Dionex solvent lines.
- 3. Place the entire assembly in the solvent reservoir.
- 4. Tighten the reservoir cap hand-tight.



Fig. 17: Connecting the solvent lines to the reservoir

When replacing a solvent line, remove the frit first, then the solvent line, and then the retaining guide.

5. The solvent supply lines are connected to the pump at the factory and routed through the opening between the front panel door and top cover.



Fig. 18: Pre-connected solvent lines on a UltiMate 3000 system pump

- ♦ All pumps (except LPG-3400 and HPG-3200P) The solvent supply lines from the reservoirs and the pump are connected to the degassing module of an appropriate SRD-3x00 Solvent Rack. For information about how to connect the tubing to the degasser, see the Solvent Rack Manual.
- ◆ *LPG-3400*

The supply lines from the solvent reservoirs and the solvent lines are connected to the degassing module of the integrated vacuum degasser.

• *HPG-3200P* The supply lines from the solvent reservoirs are connected directly to the pump inlet.

4.4 Connecting the Rear Seal Wash System

The peristaltic pump is installed at the top right in the pump. The tubing under the peristaltic pump lever remains compressed and does not relax, thus blocking the wash solution. This can happen if the pump is not running for a longer period, for example, during shipment. That is why the pump is shipped with the active rear seal wash tubing *bypassing* the peristaltic pump.

1. Verify that the peristaltic tubing is engaged in the peristaltic pump. If it is not, press the lever to the right, place the tubing in the pump, and release the lever.



Fig. 19: Peristaltic pump

- 2. Fill the liquid reservoir. Follow the steps that are appropriate for your seal wash setup.
 - a) For the standard setup of the rear seal wash system, the components are connected at the factory as shown in Fig. 7 (→ page 22). Fill the liquid reservoir as described on page 50.
 - b) If you want to perform rear seal washing from an external liquid reservoir, follow the steps on page 52.
 - **Important:** Make sure that the liquid used for rear seal washing is miscible with the solvent. This is to avoid impairing the tightness of the pump.
 - ▲ Important: Assurez-vous que le liquide utilisé pour le rinçage du joint arrière est miscible avec le solvant afin de préserver l'étanchéité de la pompe.

4.4.1 Filling the Internal Wash Liquid Reservoir

The pump is delivered with the rear seal wash system being connected as shown in Fig. 7 (\rightarrow page 22). Remove and fill the liquid reservoir:

1. Hold the liquid reservoir including the holding clip and push both parts together vertically toward the top. The holding clip disengages.



Fig. 20: Liquid reservoir of the rear seal wash system

2. While holding the reservoir by its cap, unscrew the reservoir including the holding clip from the cap.

Avoid disconnecting the tubing from the cap. If the tubing is disconnected, make sure to reconnect the tubes to the correct connection ports (\rightarrow Fig. 21). Be careful not to interchange them.



Fig. 21: Connection ports on the reservoir cap

3. Fill the reservoir. Approximately 50 mL of fluid will be sufficient. Observe the information regarding the composition of the seal-washing medium (\rightarrow page 51).

Composition of the seal-washing medium

For reliable detector performance, make sure that the seal-washing medium is slightly conductive. Standard HPLC-grade water with, for example, 5-10% isopropanol or methanol, is appropriate. If a medium other than HPLC-grade water has to be used due to the solubility of the delivered solvent, make the medium slightly conductive using the appropriate additives. (Do not use additives with a high salt content or additives that cause solid residuals upon evaporation.) Make sure that the seal-washing medium is compatible with the silicone tubing.

NP applications

For these applications, isopropanol with 0.1% sulfuric acid is recommended to ensure that the liquid is conductive.

i Tips: Regularly check the liquid level in the internal liquid reservoir, making sure that the level is always between the minimum and maximum markers on the label.

The seal-washing liquid is delivered in circles. Therefore, replace the liquid in the liquid reservoir at regular intervals.

- When using solvents *without* buffer additions, replace the liquid in the reservoir *at least* once a week.
- When using solvent *with* buffer additions or solvents with a high salt content, replace the liquid in the reservoir *at least* every second day. This is especially important with solvents with a high salt content to avoid salt concentration and prevent damage to the piston seals.

4.4.2 Installing and Filling the External Wash Liquid Reservoir

- 1. Disconnect both silicone tubes from the internal liquid reservoir (\rightarrow Fig. 7, page 22).
- 2. Fill the external liquid reservoir with an appropriate seal-washing liquid. Observe the information about the composition of the seal-washing medium (\rightarrow page 51).
- 3. Connect the silicone tubing from the peristaltic pump with the liquid reservoir. If necessary, prolong the tubing, for example, using part of the silicone tubing from the pump accessories kit.



Fig. 22: Rear seal wash system with external reservoir (here for a DGP-3600)

- 4. Place the liquid reservoir in the Solvent Rack of the UltiMate 3000 system.
- 5. Direct the tubing from the last pump block into an appropriate waste container. If necessary, prolong the tubing, for example, using part of the silicone tubing from the pump accessories kit.
- 6. Flush the system with the seal-washing liquid. Draw liquid into a syringe at the open end of the waste line.
- **I** Tips: Regularly check the liquid level in the liquid reservoir. There will be no warning when the reservoir is empty.

Replace the liquid in the liquid reservoir at regular intervals. This is especially important with solvents with a high salt content to avoid salt concentration and prevent damage to the piston seals.

4.5 Connecting the Drain Tubing

To discharge liquid leaks and waste, the pump has a drain port at the bottom right of the instrument.



Fig. 23: Drain port

Direct liquid leaks to waste through the drain system of the UltiMate 3000 system. The required components are shipped with the pump, but can be ordered also separately. The drain kit (part no. 6040.0005) includes all components for system drainage and detailed installation instructions.

4.6 Purging the Pump

Purging the pump means rinsing the system for a short time at a higher flow rate. Select one of the following alternatives:

- Purge the pump manually (\rightarrow section 4.6.1).
- If the UltiMate 3000 system includes a WPS-3000SL autosampler, you can purge the pump also via the autosampler (\rightarrow page 56).

i Tip: The two pumps of a DGP-3600 must be purged separately.

After the purge cycle has been initiated, the pump is purged with the default settings.

| Pump | Flow Rate (Purge flow) | Time (Purge time) |
|----------------------|------------------------|-------------------|
| Analytical pump | 6 mL/min | 5 min |
| Micro pump | 2 mL/min | 5 min |
| Semipreparative pump | 30 mL/min | 5 min |

You can change the default settings if necessary, either in Chromeleon or on the front panel display. In Chromeleon, change the **PurgeTime** and **PurgeFlow** settings in the **Commands** dialog box or on the control panel for the pump. On the front panel display, enter the desired values under **Purge time** and **Purge flow** on the **Preferences** menu.

4.6.1 Purging the Pump Manually

1. Attach a piece of silicone tube to the purge outlet nozzle on the outlet unit. (Silicone tubing is provided in the accessory kit for the pump.)



Fig. 24: Purge valve and purge outlet nozzle

- 2. Attach the free end of the tube to a plastic syringe. (A syringe is provided in the accessory kit for the pump.)
- 3. Loosen the purge valve by **one** turn.
- 4. To be able to generate negative pressure, an autosampler or another type of pressure drop needs to be connected to the pump outlet. If necessary, cap the pump outlet.

5. Purge the pump from Chromeleon or from the front panel display.

To purge the pump from Chromeleon

- a) In Chromeleon, open the **Commands** dialog box for the pump.
- b) Under Pump, set the channel to be purged to 100%. Chromeleon calculates the value for %A according to the following formula %A = 100% - (%B + %C + %D). Thus, %A is 100% if you set the other components of the eluent to 0%.
- c) *If the UltiMate 3000 system includes a WPS-3000SL autosampler* To avoid that the pump is inadvertently purged via the autosampler, verify that **PurgeViaSampler** is set to **No**.
- d) Start the purge cycle. To do so, set **Purge** to **On**. The purge cycle is performed with the selected settings (\rightarrow page 54).
- e) When the specified **PurgeTime** has elapsed, **Purge** is automatically reset to **Off**. You can stop purging also manually by setting **Purge** to **Off** yourself.
- f) Repeat the above steps for *all* channels channels (even if they are not used for the application) until all air bubbles are gone.

To purge the pump from the front panel display

- a) On the front panel display, select the **Control** menu.
- b) Set the channel to be purged, for example %A, to 100%
- c) Start the purge cycle. To do so, set **Purge** to **On**. The purge cycle is performed with the selected settings (\rightarrow page 54).
- d) When the specified **PurgeTime** has elapsed, **Purge** is automatically reset to **Off**. You can stop purging also manually by setting **Purge** to **Off**.
- e) Draw the liquid with the syringe.
- f) Repeat the above steps for *all* channels (even if they are not used for the application) until all air bubbles are gone.

I Tip: You can start the purge cycle also by selecting **Purge** on the navigation bar.

6. Close the purge valve. Do not overtighten the valve screw. In case of leakage, tighten more as necessary.

Do not use any tools to tighten the purge valve screw. Overtightening can destroy the cap seal. Open or close the purge valve only when the system pressure is down.

4.6.2 Purging the Pump via the WPS-3000SL Autosampler²

If the UltiMate 3000 system includes a WPS-3000SL autosampler, you can purge the pump with Chromeleon via the membrane pump of the autosampler. In this way, you can remove small amounts of air, for example, after exchanging the solvent. For an optimum performance, the purge flow should be as high as possible, for example

| Ритр Туре | Flow Rate (Purge Flow) |
|----------------------|------------------------|
| Analytical pump | 6 mL/min |
| Micro pump | 2 mL/min |
| Semipreparative pump | 30 mL/min |

In addition, the filter frit in the inline filter of the outlet unit should be permeable. If necessary, replace the filter frit (\rightarrow page 112) or purge the pump manually (\rightarrow page 54).

- 1. Open the properties of the pump in the Chromeleon Server Configuration program $(\rightarrow page 44)$.
- 2. On the **Devices** page, verify that the pump that is fluidically connected to the autosampler is selected in the **purges via WPS-3000SL** field (\rightarrow page 39).
- 3. Open the **Commands** dialog box for the pump (\rightarrow page 62) and check and change the following settings if necessary:

| Property | Setting |
|-----------------|--|
| SamplerDevice | Shows the name of the autosampler via which the pump is purged. Usually, this is the name specified under Device Name on the Devices page during the installation of the autosampler. If more than one autosampler is available in the timebase, select the autosampler to be used for pump purging from the drop-down list. |
| PurgeViaSampler | Set PurgeViaSampler = Yes to purge the pump via the autosampler selected under SamplerDevice . |

4. Start the purge cycle. To do so, set **Purge** to **On**.

This is what happens: The autosampler needle descends into the wash port, and the solvent is drawn and directed to the waste via the membrane pump of the autosampler. The purge cycle is performed with the time and flow rate settings made under **PurgeTime** and **PurgeFlow**. When the specified **PurgeTime** has elapsed, **Purge** is automatically reset to **Off**. You can stop purging also manually by setting **Purge** to **Off** yourself.

² For patent reasons, this function must not be used in the following countries: China, Germany, Great Britain, Japan, and the USA.

4.7 Equilibrating the System

Before using the pump for sample analysis, equilibrate the UltiMate 3000 system:

- 1. Pump the starting solvent through the entire system until the system is free of any other liquid composition.
- 2. Heat or cool all temperature-controlled devices, such as the column oven, to the temperature required for the application.
- 3. Set the detector wavelengths and turn on the lamps. (Wavelength calibration is performed automatically.)
- 4. Monitor the pump pressure and verify that the reading is correct for the application and is stable.
- 5. Monitor the detector signal and verify that the baseline signal is at the expected reading for your application and is stable.

Perform system equilibration in Chromeleon or select the required commands and parameters on the front panel menus of the instruments.

To equilibrate the system from Chromeleon

- Select and perform the operating commands and parameters from the **Commands** dialog box.
- Create and run an equilibration program to automate the process (\rightarrow page 65).
- Use the SmartStartup Wizard to create and run the equilibration program (see below).

To create the equilibration program with the SmartStartup Wizard

The SmartStartup Wizard in Chromeleon guides you through the process of creating and running an equilibration program.

- 1. To open the wizard, select SmartStartup on the Batch menu.
- 2. Follow the instructions as they appear on each page of the wizard. For additional information about a page, click **Help**.
- 3. After you finish the wizard, Chromeleon
 - Generates an equilibration program and sequence.
 - ◆ Opens the equilibration control panel for the instruments in the timebase (→ Fig. 25, page 58).
 - Opens the **Start Batch on** dialog box.

Click Start to begin equilibration.



The equilibration panel shows the equilibration status of each instrument in the system.

Fig. 25: Equilibration panel

To equilibrate the system from the front panel menus

Select and perform the operating commands and parameters on the front panel menus of the instruments. For information about the pump menus, see section 5.4.2 (\rightarrow page 68). For information about the menus of other system modules, see *Operating Instructions* for the respective module.

5 Operation and Maintenance

The pump can be controlled by the Chromeleon Chromatography Management System. For details, see section 5.3 (\rightarrow page 61).

In addition, function keys and menus are available on the front panel display to facilitate operation during, for example, initial installation of the pump, diagnostics, and maintenance. They allow you to perform certain actions directly from the pump display:

- Executing certain commands (purge, set flow)
- Setting parameters (leak sensor mode, brightness and contrast of the screen display)
- Viewing diagnostic information
- Viewing and changing the pump configuration

For details, see section 5.4 (\rightarrow page 67).

5.1 Power-Up

To start the pump for the first time, turn on the main power switch on the rear panel of the pump. The following sequence of events occurs when the pump is powered up:

- For a short time, general information about the pump appears on the front panel display: device type, serial number, bootloader and firmware versions.
- The pump runs a series of internal tests. During these self-diagnostics, all of the main components are checked. When testing is complete and has been successful, the initial screen changes to the status screen (→ page 60).
- If an error is detected, the pump is not ready for analysis. The **Status** LED on the front panel door changes to red and a message appears on the front panel display. If the pump is operated with Chromeleon, the message is logged also in the Chromeleon Audit Trail. Turn off the pump, take appropriate remedial action (→ page 90), and turn on the pump again.

For routine operation, leave the main power switch on. For routine on/off control, use the standby button on the front of the pump (\rightarrow page 17). Press and hold the button for one second to allow the pump to change the mode. Turn the main power switch off when instructed to do so, for example, before performing a service procedure.

5.2 Status Screens

When the self test was successful, the initial screen changes to the status screen. The pump type determines the appearance of the screen:



Fig. 26: Example of a status screen

The status screen shows the following information:

• Pump name

The pump name is the name specified on the **Devices** page in Chromeleon Server Configuration program (\rightarrow page 39).

- Flow
- Pressure
- Components of the solvent in percent of the total flow

In addition, the following information may appear on the screen:

| The screen shows | When |
|-------------------------------|---|
| On or Flow on | The pump is delivering with the specified flow rate. |
| Off or Flow off | The pump is not delivering. |
| Purge | The pump is purged. |
| Master flow or Column flow | The UltiMate 3000 system includes a DGP-3600M(B), LPG-3400M(B), or HPG-3x00 pump and an FLM-3x00 Flow Manager. The values refer to the flow rate and pressure before the flow splitter (Master flow) or after the flow splitter (Column flow). For more information, see page 77. |

You can adapt the screen brightness and contrast to your requirements if necessary (\rightarrow page 85).

5.3 Operation with Chromeleon

Before you begin, verify that

- 1. The Chromeleon software is installed on the computer and the license code is entered.
- 2. The pump is connected to the Chromeleon computer via a USB connection.
 - **i** Tip: Verify that Chromeleon is installed on the computer and that the license code is entered *before* you connect the pump to the USB port on the Chromeleon computer and turn on the pump power. Only then, the USB driver for the pump is automatically loaded and the Windows operating system can detect the pump when the power is turned on.
- 3. The pump is set up in Chromeleon, as described in section 3.5 (\rightarrow page 34).

Before you can operate the pump with Chromeleon, you have to connect the timebase in which the pump is installed to the Chromeleon client program (\rightarrow page 61).

Two modes of software control are available:

- Direct control with the parameters and commands in the **Commands** dialog box (F8 box) or on a control panel (→ page 62)
- Automated control with a control program (PGM) (\rightarrow page 65)

5.3.1 Connecting to Chromeleon

- 1. Start the Chromeleon Server Monitor if it is not already running (\rightarrow page 34).
- 2. Start the Chromeleon client by clicking the Chromeleon icon $\overset{\bullet}{\Longrightarrow}$ on the desktop.

If the Chromeleon icon is not on the desktop, click **Start** on the taskbar, point to **Programs** (or **All Programs**, depending on the operating system), point to **Chromeleon**, and then click **Chromeleon**.

3. Connect the Chromeleon client to the timebase in which the pump is installed. For details about how to do this from the **Commands** dialog box, see page 62. For details about how to do this on a control panel, see page 64.

When the pump is correctly connected to Chromeleon:

- The **Connected** LED on the front panel is green.
- Front panel input is disabled.
- Functions for estimating the lifetime of consumables and monitoring and recording service and (re)qualification information are provided (\rightarrow page 86).

- Diagnostics functions are available for checking the performance of certain pump components as well as the overall performance of the instrument (\rightarrow page 87).
- The **Standby** button on the front panel remains active.

Before turning off the pump by the main power switch, always **disconnect** the module in Chromeleon.

5.3.2 Direct Control

With direct control, you select operating parameters and commands in the **Commands** (F8) dialog box. Direct commands are executed as soon as they are entered. For routine operation, most parameters and commands are available also on a control panel.

To open the Commands dialog box for the pump

- 1. Open a control panel (any panel is possible). To open a control panel, open the Chromeleon Browser and double-click a control panel in the **Dionex Templates/Panels** folder.
- 2. Connect the control panel to the timebase in which the pump is installed. On the **Control** menu, select **Connect to Timebase**, and then select the timebase on the **Timebase** tab. For information about the **Timebase** dialog, click **Help**.

Tip: The **Control** menu is visible only when a control panel is already open.

- 3. Press the F8 key or select **Command** on the **Control** menu.
- 4. To see the parameters and commands that are available for the pump, click the plus sign next to **Pump**.
 - **Tip:** For a DGP-3600, separate entries for the right and left pump appear under the main **Pump** entry. The pump version (A or M) determines which entries appear. These are **PumpRight** or **MicroPump** for the right pump and **PumpLeft** or **LoadingPump** for the left pump.
The commands and parameters available in the dialog box vary, depending on the

- Chromeleon version
- options selected in the **Properties** dialog for the pump (\rightarrow page 37).
- display filter level (Normal, Advanced, or Expert)
- 5. Change the display filter level if necessary. Right-click in the commands list and select the filter level on the menu.



Fig. 27: Commands dialog box

6. Verify that the pump is connected to Chromeleon. If it is not, select **Connect** to connect the pump.

For a list of the commands and properties that are supported for the pump, see the Chromeleon Help. In addition to the pump commands and parameters, the **Commands** dialog box provides access to all of the commands and parameters available for all devices that are installed in the selected timebase.

Tip: If the UltiMate 3000 system includes an FLM-3x00 Flow Manager and if you specified on the **Configuration** page for the flow manager that the column pressure is recorded as a separate channel, a **ColumnPressure** entry appears in the **Commands** dialog box (under the related **ColumnOven**) in the addition to the pump pressure entries (\rightarrow page 40).

To open a control panel

1. On the **View** menu, click **Default Panel Tabset** or click the corresponding icon on the toolbar **10**, and then connect to the Chromeleon server.

Chromeleon creates centralized control panels, called panel tabsets (\rightarrow Fig. 28, page 64), for all timebases available on the Chromeleon server. A panel tabset provides control panels for the individual instruments in a timbase and, in addition, one or more panels for performing system-wide functions, for example, creating and running sequences. For more information about panel tabsets, see the Chromeleon Help.

- 2. On the **Panel Tabset** for your timebase, click the **Pump** page (\rightarrow Fig. 28).
- 3. Verify that the pump is connected to Chromeleon (the LED next to the **Connect** button is green). If it is not, click **Connect**.



Fig. 28: Pump control panel on the panel tabset

The control panel provides access to the operating parameters and commands required for routine operation of the pump. Additional functions are available in the **Commands** dialog box. To open the **Commands** box from the panel tabset, select **Command** on the **Control** menu.

5.3.3 Automated Control

With automated control, you create a program file (PGM) for automated operation of the pump. Programs can be created automatically with the help of a software wizard or manually by editing an existing program.

In addition to programs for sample analysis, you can also create programs for special purposes, for example, to automate system shutdown (\rightarrow page 91) or to ensure that the system automatically restarts operation as desired after a power failure. For details, see the Chromeleon Help.

To create a program with the Program Wizard

1. On the File menu, select New, and then select Program File.

The wizard guides you through program creation. On each wizard page, make the desired settings or accept the default values. For additional information about a page, click **Help**.

- 2. After you finish the wizard, Chromeleon automatically creates the corresponding program.
- 3. To start the program, follow the steps on page 66.

To create a program manually

1. Open an existing program.

Select and double-click the program you want to open.

—Or—

On the **File** menu, select **Open**. In the dialog box, select **Program** on the **Object of Type** list and select the program.



Fig. 29: Chromeleon program file (here: program shown in the Commands view)

2. Change the settings in the program as desired.

The easiest way is to edit a program is to do this in the Device Views (\rightarrow Fig. 29). Click a device icon and change the settings on the device pages. Editing the program in the Device Views ensures correct command syntax.

If you cannot edit a certain parameter in the Device View, click **Commands** to open the Commands View. The Commands View shows the entire program, listing the control commands in chronological order. Edit the parameter of interest or enter a new parameter. For more information, see the Chromeleon Help.

3. To start the program, follow the steps below.

To start a program

Program for sample analysis

- 1. Create a sample list (sequence). A sequence must include the program and a method for evaluating the sample data (for example, for peak identification, area determination, and amount determination).
- 2. Assign the program and method to each sample on the list.
- 3. Add the sequence to the batch and start the batch.

For information about each of the above steps, see the Chromeleon Help.

Other programs

Add the program to the batch and start the batch.

5.4 Display Screens (Function Keys and Menus)

Via the function keys and menus that are available on the front panel display, you can make several settings or execute certain commands directly from the pump. For information about the function keys, see section 67 and page 69. For information about the menus, see section 5.4.2 (\rightarrow page 68).

5.4.1 Showing the Function Keys

Four white spots on the front panel mark the positions of four function keys: **Menu**, **Flow on**, **Set flow**, and **Purge** (\rightarrow table on page 67). To show the keys, select the position of the utmost left spot on the front panel display with the menu pen (part no. 6300.0100). The menu pen is included in the pump shipment.



Fig. 30: Showing the function keys

The function keys replace the information in the bottom line of the status screen. If no key is selected, the bottom line of the status screen is restored after about 5 seconds.



Fig. 31: Function keys

| То | Select |
|--|----------|
| Open the Main menu (\rightarrow page 70). | Menu |
| Have the pump deliver with the specified flow rate. The name of the key changes to Flow off when the pump is delivering with the specified flow rate. Select Flow off to have the pump stop delivering. | Flow on |
| Enter the pump flow. | Set flow |
| Purge the pump. The channel set on the Control menu (\rightarrow page 71) is purged with the Purge flow and Purge time values from the Preferences menu (\rightarrow page 72) being considered. For more information about purging the pump, see section 4.6 (\rightarrow page 54). | Purge |

Front panel input is disabled while the pump is connected in Chromeleon.

5.4.2 Pump Menus

Fig. 32 shows an overview of the pump menus. For information about the general menu layout and structure, see page 69. For information about the commands and parameters that are supported by the menus, see sections 5.4.2.2 through 5.4.2.6 (\rightarrow page 70 and following pages).



Fig. 32: Menu structure

5.4.2.1 General Menu Layout and Structure

In general, the menu layout is as follows:



Fig. 33: Menu layout (here: Control menu)

| No. | Description |
|-----|---|
| 1 | Reports the menu name and the number of items on the menu list. |
| 2 | The menu items appear on a list and are numbered consecutively. The selected item is underlined. |
| 3 | Navigation bar |

Select an item with the arrow up or down key—the selected item is underlined. Confirm your selection with **Select**. **Back** returns you to the previous menu level.

The selected menu item or parameter determines which keys appear on the navigation bar:

| То | Select |
|---|--------|
| Return to the previous entry on a list. | ^ |
| list, after reaching the first line (\rightarrow Key autorepeat, page 74). | |
| Increment numerical values. | ^ |
| Proceed to the next entry on a list. | \vee |
| If the list contains 5 or more items, you can use the arrow down key to scroll down through the list, after reaching the forth line (\rightarrow Key autorepeat, page 74). | |
| Decrement numerical values. | \vee |
| Proceed to the next figure in a number. Any decimal point is skipped. | > |
| Confirm the selection and activate the input field if applicable. If an item is read-only, the Select key will not be available. | Select |
| Return to the previous menu level. | Back |
| Confirm the selection or input. | OK |
| Cancel the action and restore the last value. | Cancel |
| Note: Depending on the selected option, specific keys may replace these general keys. | |

If an error is found, one or more messages appear on the front panel display. In this case, the **Prev**, **Next**, and **Clear** keys appear on the navigation bar.

| То | Select |
|------------------------------------|--------|
| Return to the previous message. | Prev |
| Proceed to the next message. | Next |
| Remove a message from the display. | Clear |

5.4.2.2 Main Menu

The Main menu provides top-level access to the menu structure. To open the Main menu, show the function keys and select Menu (\rightarrow page 67).



Fig. 34: Main menu

For details about the menus, see the following sections:

- Control Menu (\rightarrow page 71)
- Preferences Menu (\rightarrow page 71)
- Diagnostics Menu (\rightarrow page 73)
- Configuration Menu (\rightarrow page 74)

5.4.2.3 Control Menu

On the **Control** menu, you can set the flow rate, the components of the solvent, and the pressure limits. You can start or stop the pump flow and purge the pump. In addition, you can move the pump into the appropriate position for piston installation.

| > Control (112) | | |
|-----------------------------|--------|------|
| <u>1. Flow</u> | | |
| 2. Flow rate | | |
| 3. %A | | |
| 4. %B | | |
| 5. %C | | |
| 6. %D | | |
| 7. Min pressure | | |
| 8. Max pressure | | |
| 9. Selected pump | | |
| 10. Purge | | |
| 11. Change right pump pisto | ns | |
| 12. Change left pump piston | S | |
| | Select | Back |

Fig. 35: Control menu

| То | Select |
|--|---|
| Start the pump flow with the selected flow rate or to stop the flow. | Flow |
| Set the flow rate. | Flow rate |
| Enter the related component of the solvent in percent of the total flow. | %A through %D |
| Set the maximum pressure. | Max. pressure |
| Set the minimum pressure. | Min. pressure |
| Specify for which pump of a DGP-3600, the values can be entered or displayed and the commands are executed. | Selected pump |
| Purge the pump. The selected channel is purged with the Purge flow and Purge time values from the Preferences menu (\rightarrow page 72). For more information about purging the pump, see section 4.6 (\rightarrow page 54). | Purge |
| Move the related pump into the appropriate position for piston installation $(\rightarrow page 127)$. | Change right pump pistons Change left pump pistons |

5.4.2.4 Preferences Menu

On the **Preferences** menu, you can make the basic settings for the pump.



Fig. 36: Preferences menu

| То | Select |
|--|-------------------|
| Turn the rear seal wash system on (Interval or Automatic) or Off . Interval activates rear seal washing once per rear seal wash interval for five minutes. However, the detector is <i>not</i> active. This means that leakage monitoring of the piston seals is disabled. Automatic periodically activates rear seal washing once per rear seal wash interval until the detector has counted 30 drops. The detector is active, that is, piston seal leakage is monitored. <i>Tip:</i> The default rear seal wash interval is 60 minutes. If the pump is operated with Chromeleon, you can change this setting via RearSealWashInterval (\rightarrow page 81). | Rear seal wash |
| Specify the leak detection threshold for the rear seal wash system. If the limit is met, a message appears on the front panel display. | Rear seal limit |
| Turn the degasser of an LPG-3400 or SRD-3x00 Solvent Rack on or off. | Degasser |
| Set how long the pump is purged. The default purge time is 300 seconds. | Purge time |
| Set the flow rate for purging. The allowed range depends on the pump type. The default purge flow is as follows: 6 mL/min for analytical pumps 2 mL/min for micro pumps 30 mL/min for semipreparative pumps. | Purge flow |
| Set the upper limit for the flow rate acceleration. For more information, see page 78. | Flow acceleration |
| Set the upper limit for the flow rate deceleration. For more information, see page 78. | Flow deceleration |

5.4.2.5 Diagnostics Menu

The **Diagnostics** menu provides information for diagnostic purposes (read-only). In addition, you can perform a self test.



Fig. 37: Diagnostics menu

| То | Select |
|--|---|
| Perform a self test. If an error or mechanical fault is detected, the Status LED on the front panel changes to red and a message appears on the font panel display. | Self test |
| See the compression value of the last stroke. Note: If the value is very high and the pressure is low, the pump might not be purged completely. | Compression right <i>or</i> Compression left |
| See the workload of the related pump block or pump of a DGP-3600 since the pump has been operated for the first time. The workload is calculated from the flow rate, pressure, and time. | Workload right or Workload left |
| See the pump model. | Model |
| See which firmware version is installed. | Firmware version |
| See the bootloader version. | Bootloader version |
| See the serial number of the pump. | Serial number |
| See the operating hours of the pump since the pump has been turned on for the first time. | Total operating hours |
| See the operating hours of the pump since the pump has been turned on. | Actual operating hours |

5.4.2.6 Configuration Menu

The **Configuration** menu provides information about the pump configuration and allows you to make the required settings and change the settings:



Fig. 38: Configuration menu

| То | Select |
|--|---------------------------|
| Set the display and function key parameters Brightness —Sets the screen brightness. The input is in per cent. Contrast —Sets the screen contrast. The input is in per cent. Key sound —Sets whether an acoustic beep sounds when you select a function key: On —yes or Off —no. Key autorepeat — sets whether the arrow up and arrow down keys can be used to scroll through the items on a menu list if the list contains 5 or more entries: On —yes or Off —no. | Display & soft keys |
| Set whether leak detection is be performed and how you are alerted in case of an alarm: Enabled —activates leak detection. When a leak is detected, a message appears on the front panel display and an acoustic beep sounds. Silent —activates leak detection. When a leak is detected, a message appears on the front panel display, but <i>no</i> beep sounds. Disabled —deactivates leak detection | Leak sensor mode |
| Set the pressure unit. | Pressure unit |
| See the system pressure (\rightarrow page 20). For a DGP-3600 pump, the system pressure is the pressure of the right pump. | DAC output |
| See whether the values on the status screen refer to the flow rate and pressure <i>before</i> the flow splitter (Master flow) or <i>after</i> the flow splitter (column flow) when the UltiMate 3000 system includes an DGP-3600M(B), LPG-3400M(B), or HPG-3x00A or M, and an FLM-3x00 Flow Manager. | Display mode |
| Restore the factory settings. In the Reset to factory defaults? dialog box, select OK to confirm the restore or select Cancel to keep your settings. | Reset to factory defaults |

5.5 Operational Settings

5.5.1 Choosing the Solvents

All parts of the pump that may be exposed to solvents are made, for example, of stainless steel, titanium, PCTFE, PTFE, PEEK, or sapphire (\rightarrow page 159). For information about the chemical resistance of PEEK, see the table in section 12.1 (\rightarrow page 175).

• Use only standard solvents and water of at least HPLC grade or better LC-MS grade (0.2 µm, filtered), and buffers that are compatible with the flow path materials.

If water from water purification systems that are not property maintained is used, polymeric contamination may seriously damage the column and rapidly block solvent frits.

- Mind the special properties of the solvents, such as viscosity, boiling point, UV absorption (UV/VIS detector), refractive index (refractive index detector), and dissolved gas (degasser).
- Stainless steel pumps

Buffer concentration: Typically up to 1 mol/L (< 0.1 mol/L chloride ions). Make sure to use the active rear seal wash system (\rightarrow page 49).

pH range: 1 through 13.

When using filter frits with a porosity of 0.5 μ m, Thermo Fisher Scientific recommends installing PEEK filter frits for pH values < 4.

- Make sure to use special (highly pure) solvents. They are usually labeled accordingly by the vendor.
- **Important:** Thermo Fisher Scientific advises against recycling the solvents or using methanol from aluminum reservoirs. This may impair the performance of the seals.
- ▲ Important: Thermo Fisher Scientific déconseille de recycler les solvants ou employer du méthanol stocké dans des réservoirs en aluminium. Ceci peut affecter les performances des joints.

The pump is shipped with reversed phase piston seals. You can replace these seals with normal phase seals if required. For information about the installation procedure and the corresponding part numbers, see section 7.5.2 (\rightarrow page 119).

5.5.2 Linking the Pump to the Autosampler

Only if the UltiMate 3000 system includes an UltiMate 3000 series autosampler

If the UltiMate 3000 system includes an UltiMate 3000 series pump and a WPS-3000PL, WPS-3000SL, or ACC-3000 autosampler, you can specify to which pump the autosampler is linked. Thermo Fisher Scientific recommends always specifying a pump because this setting

- Indicates which pump delivers the flow. The information is important if you want to operate the system in bypass mode (possible only with WPS-3000SL). For information about the bypass mode, see the *Autosampler Manual* or Chromeleon Help.
- Allows synchronizing the injection command of the autosampler with the strokes of a low-pressure gradient pump (DGP-3600 or LPG-3400). Synchronization ensures that all injections are performed at the same phase of the pump cycle, and thus help enhancing considerably the retention time precision with gradient applications.
- 1. Start the Server Configuration program (\rightarrow page 36).
- 2. Right-click the autosampler in the timebase and select **Properties** on the menu.
- 3. On the **Segments / Pump Link** page, select the pump to which the autosampler is linked on **Flow through sampler is delivered by pump(s)** list.
 - ♦ If the UltiMate 3000 system includes a DGP-3600 pump Select UM3PUMP_L_STRK to link the autosampler to the left pump. To link the autosampler to the right pump, select UM3PUMP_R_STRK.
 - If the UltiMate 3000 system includes a pump other than a DGP-3600 Select UM3PUMP_STROKE.
 - If you do not want to link the autosampler to a pump Select <None>.

The pump link setting on the **Segments / Pump Link** page is the default setting. To change the setting for a specific application, use the **SyncWithPump** and **PumpDevice** properties in the **Commands** dialog box or in the program (PGM).

- To disable synchronization, set SyncWithPump to Off.
- To use a different pump, enter the pump under **PumpDevice**, and then verify that **SyncWithPump** is set to **On**.

The default setting on the Segments / Pump Link page remains unchanged.

5.5.3 Linking the Pump to the Flow Manager and Setting the Display Mode

Only if the UltiMate 3000 system includes a micro pump or high-pressure gradient pump and an FLM-3x00 Flow Manager

If the UltiMate 3000 system includes one of the following pumps

- DGP-3600M or DGP-3600MB
- LPG-3400M or LPG-3400MB
- HPG-3200A or HPG-3200M
- HPG-3400A oder HPG-3400M

and an FLM-3x00 Flow Manager, you can connect the pump with the flow splitter in the Flow Manager.

When you assign a flow splitter to a pump, the pump flow settings apply directly to the flow on the splitter outlet, that is, the split ratio is considered automatically. In the standard configuration of the UltiMate 3000 system with a DGP-3600M or DGP-3600MB, Thermo Fisher Scientific recommends assigning the flow splitter to the right pump (MicroPump).

- 1. Start the **Server Configuration** program (\rightarrow page 36).
- 2. Right-click the pump in the timebase and select **Properties** on the menu.
- 3. On the **Devices** page, assign the flow splitter in the uses FLM-3x00 Flow Splitter field (\rightarrow page 41).
- 4. Specify which information is displayed on the front panel display of the pump: Flow rate and pressure before the flow splitter (= master flow) or flow rate and pressure after the flow splitter (= column flow).

Open the **Commands** dialog box for the pump. Click **DisplayMode** and select **Master** or **Column** from the list. As an alternative, you may set the display mode on the **Configuration** menu on the pump display.

5.5.4 Setting the Flow Rate, Flow Acceleration, and Flow Deceleration

In Chromeleon or on the pump menus, you can set how fast the pump starts delivering with the selected flow rate (flow acceleration) and how fast the pump flow is reduced (flow deceleration).

- If the values are too low, it will take accordingly long for the pump to reach the necessary flow, and thus to build up the necessary pressure or reduce the flow and thus the pressure as required.
- If the values are too high, this may destroy the filter frit in the inline filter of the pump or reduce the lifetime of the column.

Thermo Fisher Scientific recommends that you set these parameters to values between 1/3 and factor 3 of the (column) flow rate. If you set the parameters to values outside this range, a warning appears in Chromeleon.

To set the flow rate, flow acceleration, and flow deceleration from Chromeleon

- 1. Open the **Commands** dialog box for the pump.
- 2. Select **Flow** and enter the desired flow rate.

The allowed flow rate range is indicated in the properties of the pump, depending on the pump type on the **Head Type & Limits** page or on the **Right Pump** and **Left Pump** pages (\rightarrow page 42). You can change the upper and lower limits for the flow rate within the allowed range and enable the **Double flow** mode if the pump is a high-pressure gradient pump (\rightarrow page 89).

3. Under **MaximumFlowRampUp**, check and change the flow acceleration setting if necessary.

Under **MaximumFlowRampDown**, check and change the flow deceleration setting if necessary.

To set the flow rate, flow acceleration and flow deceleration from the front panel display

- 1. On the front panel display, select the **Control** menu.
- 2. Select **Flow rate** and enter the desired value.

I Tip: You can enter the flow rate also by selecting **Set flow** on the navigation bar.

- 3. On the front panel display, select the **Preferences** menu.
- 4. Check and change the settings for **Flow acceleration** and **Flow deceleration** if necessary.

5.5.5 Setting the Pressure Limits

The pump firmware and Chromeleon provide default values for the upper and lower pressure limits. The limits depend on the pump type. You are free to change the limits within the allowed pressure range.

If the pump pressure is outside the specified limits, the related message appears on the front panel display. If the pump is operated with Chromeleon, the message appears also in the Chromeleon Audit Trail. In addition, Chromeleon stops the flow and aborts the batch. Check the Troubleshooting section for a short description of possible causes along with recommended courses of action (\rightarrow page 90).

• Min. pressure/Lower Limit

Helps to prevent the pump, and thus the column, from running dry. A typical setting is 10 bar.

• Max. pressure/Upper Limit

Helps to protect the column for too high a pressure. The application and column type determine the setting.

To set the pressure limits in Chromeleon

A To set the default limits

The allowed pressure range is indicated in the properties of the pump, depending on the pump type on the **Head Type & Limits** page or on the **Right Pump** and **Left Pump** pages (\rightarrow page 42). You can change the upper and lower pressure limits within the allowed range and select the pressure unit.

Tip: Changing the settings for a specific application in the **Commands** dialog box, in a program file (PGM), or on a control panel, will not change the default settings on the configuration pages.

B To change the pressure limits for a specific application

- 1. Open the **Commands** dialog box for the pump.
- 2. Select **Pressure** and enter the new limits under **LowerLimit** and **UpperLimit**.

To set the pressure limits on the front panel display

- 1. On the front panel display, select the **Control** menu.
- 2. Check and change the settings for Max. pressure and Min. pressure if necessary.

You can change the **Pressure unit** on the **Configuration** menu.

5.5.6 Recording the Pump Pressure

On the **Devices** page (\rightarrow page 39), the **Pressure Signal(s)** check box (\rightarrow page 40) is selected by default when the pump is installed and configured in Chromeleon. With this setting, Chromeleon generates the appropriate channels for recording the pump pressure. Thermo Fisher Scientific recommends recording the pump pressure in the Chromeleon program. If a problem occurs, the pump pressure channel can provide helpful information to identify and eliminate the source for the problem.

The channel is available in the **Commands** dialog box for the pump.

For all pumps (except the DGP-3600), the channel name is generated from the entry in the **Device Name** input field on the **Devices** page and is extended by **_Pressure**. For the DGP-3600, the names are generated from the entries in the **Right Pump** and **Left Pump** input fields and are extended by **_Pressure**.

In all cases, the recorded pressure is the column pressure.

Tip: If the LPG-3400M(B) or DGP-3600M(B) micro pump has the flow splitter of an FLM-3x00 Flow Manager assigned, the name ends in **_MasterPressure**. This channel records the *pump pressure* (= the pressure *before* the flow splitter).

To record the *column pressure* (= pressure *after* the flow splitter), verify on the on the Components page of the Flow Manager properties that the Pressure Signal check box is selected. The associated **ColumnPressure** channel is then available in the **Commands** dialog box under **ColumnOven**.

5.5.7 Activating Rear Seal Washing

- ▲ **Important:** If you use solvents with a high salt content, do not operate the pump without rear seal washing for a longer time (> 5 minutes). This may cause damage to the piston seals and the piston.
- ▲ Important: Si vous utilisez des phases mobiles avec une forte teneur en sel, ne faites pas fonctionner la pompe sans rinçage du joint arrière pendant un temps prolongé (> 5 minutes). Ceci peut endommager les joints de piston et le piston.

The configuration of the seal wash system determines the settings to be used and what happens:

- For the standard setup (\rightarrow Fig. 7, page 22), see the information below.
- For the alternative setup (\rightarrow Fig. 8, page 22), see page 83 for details.

For information about the seal-washing medium and about how to fill the liquid reservoir, see section 4.4 (\rightarrow page 49).

Standard setup

To activate seal washing from Chromeleon

- 1. In Chromeleon, open the **Commands** dialog box for the pump.
- 2. Click **RearSealWashSystem** and select **Automatic** on the list. Seal washing is periodically activated once per **RearSealWashInterval** until the detector in the reservoir cap has counted 30 drops.
- 3. Under **RearSealWashInterval**, check and change the time interval after which seal washing is activated. The default setting is 60 minutes.
- 4. Under **RearSealLeakage.Limit**, check and change the value for the leak detection threshold if necessary.

To activate seal washing from the front panel display

- 1. On the front panel display, select the **Preferences** menu.
- Set Rear seal wash to Automatic. Seal washing is periodically activated once per rear seal wash interval until the detector in the reservoir cap has counted 30 drops. The default rear seal wash interval is 60 minutes. If the pump is operated with Chromeleon, you can change the time interval via RearSealWashInterval.
- 3. Under **Rear seal limit**, check and change the value for the leak detection threshold if necessary.

This is what happens ...

When rear seal washing is activated, the detector in the cap on the liquid reservoir (drop counter) is active, monitoring the performance of the seal wash system, and thus detecting possible leakage of the main piston seal.

Rear seal washing performs correctly

During the delivery period of the peristaltic pump, liquid reaches the detector on the liquid reservoir (the drops are counted). This means that the seal wash system performs correctly. The liquid reservoir contains enough liquid for seal washing, the tubing is all right, and the peristaltic pump works correctly.

Malfunctioning of the rear seal wash system

If no drops reach the detector on the liquid reservoir after maximum five minutes although the peristaltic pump is pumping, this indicates that either the liquid reservoir is empty or the peristaltic tubing is broken or sticks together.

This malfunctioning may also occur if the detector on the liquid reservoir is very dirty. The following message appears "The rear seal leak sensor is malfunctioning." Clean the electrodes of the detector, using water or solvent. Be careful not to bend the electrodes.

In all cases, the following message appears in the Chromeleon Audit Trail: "Rear seal wash system has run out of wash solution". Refill the liquid reservoir or replace the defective peristaltic tubing. For more information regarding this message, see the Troubleshooting section (\rightarrow page 90).

Possible leakage of the main piston seal

If, during the period in which the peristaltic pump is idle, more drops are counted than specified under **Rear seal limit** on the **Preferences** menu (\rightarrow page 72) or under **RearSeal Leakage.Limit** in Chromeleon, this indicates possible leakage of the main piston seal. The following message appears in the Chromeleon Audit Trail "The rear seal leak count is xx (= number of counted drops) and has exceeded the limit of yy (= leak threshold value)".

Take the following remedial action:

- Visually inspect the pump for liquid leaks from the piston seals (\rightarrow page 117) or run the related leak test in Chromeleon (\rightarrow page 87).
- Replace the piston seals and support rings as necessary (\rightarrow page 123)
- Increase the value for the rear seal limit. However, increase the value only if the leakage does not affect the accuracy or precision of the retention times.

- **Tip:** The seal-washing liquid is delivered in circles. Therefore, replace the liquid in the liquid reservoir at regular intervals:
 - When using solvents without buffer additions or solvents with a high salt content, replace the liquid in the reservoir *at least* once a week. This is especially important with solvents with a high salt content to avoid salt concentration and prevent damage to the piston seals.
 - When using solvent with buffer additions, replace the liquid in the reservoir *at least* every second day.

Alternative Setup

To activate seal washing from Chromeleon

- 1. In Chromeleon, open the **Commands** dialog box for the pump.
- Click RearSealWashSystem and select Interval on the list. Seal washing is periodically activated once per RearSealWashInterval for five minutes.
- 3. Under **RearSealWashInterval**, check and change the time interval after which seal washing is activated. The default setting is 60 minutes.

To activate seal washing from the front panel display

- 1. On the front panel display, select the **Preferences** menu.
- 2. Set **Rear seal wash** to **Interval**. Seal washing is periodically activated once per rear seal wash interval for five minutes. The default rear seal wash interval is 60 minutes. If the pump is operated with Chromeleon, you can change the time interval via **RearSealWashInterval**.

This is what happens ...

As the internal liquid reservoir of the pump is not connected in this setup, the drop detector of the seal wash system is not active either. Thus, possible leakage of the main piston seal will not be monitored.

Monitor the retention times to identify possible leakage at an early stage. Stable retention times indicate that the main piston seals seal sufficiently tight. Bad retention time precision and accuracy may be due to liquid leaks from the main piston seals.

In this case, take the following remedial actions:

- Visually inspect the pump for liquid leaks from the main piston seals (\rightarrow page 117) or run the related leak test in Chromeleon (\rightarrow page 87).
- Replace the piston seals and support rings if necessary (\rightarrow page 123).

Tips: Regularly check the liquid level in the external liquid reservoir. There will be no warning when the reservoir is empty.

In this mode, the seal-washing liquid is not delivered in circles. However, replace the liquid in the liquid reservoir at regular intervals nevertheless (*at least* once a week). This is especially important with solvents with a high salt content to avoid salt concentration and prevent damage to the piston seals.

5.5.8 Purging the Pump

Select one of the following alternatives to remove air bubbles from the system:

- Purge the pump manually (\rightarrow page 54).
- If the UltiMate 3000 system includes a WPS-3000SL autosampler, you can purge the pump also via the autosampler (\rightarrow page 56).

Tip: To remove all air bubbles from the system, all channels of the pump must be filled with liquid even if one or more channels are not used for the application.

5.5.9 Activating Leak Detection

Only with the standard setup of the rear seal wash system

You can activate and deactivate leak detection in Chromeleon and on the front panel display. The running analysis will be aborted when the leak sensor reports a leak.

To activate leak detection in Chromeleon

- 1. In Chromeleon, open the **Commands** dialog box for the pump.
- 2. Select one of the following alternatives:
 - a) To activate leak detection, select LeakSensorMode, and then select Enabled or Silent. When the leak sensor reports a leak, the Status LED on the front panel changes to red. In addition, when the setting is
 - Enabled, a message appears on the front panel display and a beep sounds.
 - Silent, a message appears on the front panel display but no beep sounds.
 - b) To deactivate leak detection, select **Disabled**.

To activate leak detection on the front panel display

- 1. Show the function keys and select **Menu** (\rightarrow page 60).
- 2. Select the **Configuration** menu and select **Leak sensor mode**.
- 3. Select Enabled, Silent, or Disabled. (For details about these settings, see above.)

5.5.10 Adjusting the Screen Brightness or Contrast

You can adjust the screen brightness or screen contrast to your requirements from Chromeleon or on the front panel display.

To adjust the settings from Chromeleon

- 1. In Chromeleon, open the **Commands** dialog box for the pump.
- 2. Select **Brightness** and change the value for the screen brightness as appropriate. Select **Contrast** and change the value for the screen contrast as appropriate.

To adjust the settings on the front panel display

- 1. Show the function keys and select Menu.
- 2. Select the **Configuration** menu and select **Display & soft keys**.
- 3. Select **Brightness** and change the value for the screen brightness as appropriate. Select **Contrast** and change the value for the screen contrast as appropriate.

5.5.11 SmartStartup and SmartShutdown

The **SmartStartup** wizard (\rightarrow page 57) assists you in automating regular routine tasks. With SmartStartup, the different modules of the UltiMate 3000 system are turned on automatically and in a controlled manner. For example, SmartStartup can purge the pump of the HPLC system automatically, flush the column, and perform system equilibration. Important module parameters, such as the pressure pulsation of the pump, are monitored. When the modules operate within the specified limits, the sample sequence, which was set up before, can be started automatically. SmartStartup can be used at any time.

If you have to interrupt system operation, use the **SmartShutdown** wizard to create a program to set the HPLC system into standby mode or to automate shutdown of the system $(\rightarrow page 91)$.

5.6 Special Chromeleon Functions

This section provides a short overview of some special functions that Chromeleon supports for the pump. All of these functions are available in the **Commands** dialog box (unless otherwise noted). In addition, some functions are available also on the control panel for the pump. For additional information about a function, see the Chromeleon Help.

5.6.1 Predictive Performance

Predictive Performance provides various functions for estimating the lifetime of consumables and for monitoring and recording service and (re)qualification information.

Commands Dialog Box

Open the **Commands** dialog box for the pump and enter the limits for the predictive performance parameters. For a complete list of available commands and parameters, see the Chromeleon Help. To keep the predictive performance information up-to-date, Thermo Fisher Scientific recommends the following actions:

| After you have | Perform the following command |
|--|-------------------------------|
| Replaced the check valve | CheckValveServiceDone |
| Replaced the filter frit in the inline filter of the outlet unit | OutletFritChanged |
| Replaced the piston | PistonsChanged |
| Replaced the piston seal | SealsChanged |
| Replaced the peristaltic pump tubing | RearSealWashTubeChanged |
| Serviced the pump (for example, annual maintenance) | ServiceDone |
| Performed instrument qualification | QualificationDone |

These commands reset the related counters and update the information when the action was performed.

Control Panel

On the control panel for the pump, click **Wellness**, **Qualification**, and **Service** to see the related predictive performance commands and parameters on separate panels. On these panels, you can enter the limits and reset the counters. In addition, wellness bars provide visual indicators of qualification and service periods (Chromeleon 6.80 and later). The color-coding of the wellness bars provides information about the status:

| Color | Description |
|--------|--|
| Green | OK. |
| Yellow | The value will soon reach the specified limit and/or the related component needs servicing or should be replaced soon. |

| Color | Description |
|--------|---|
| Orange | (Only for monitoring Qualification properties) The value has reached the specified limit. However, a Grace Period has been specified during which the component may still be operated. |
| Red | The value has reached the specified limit or the specified grace period has expired. Replacement, servicing, or qualification of the component is overdue. The component can no longer be operated and it is not possible to start a batch. |

In addition, a message appears in the Chromeleon Audit Trail when a limit has been reached.

5.6.2 Pump Diagnostics

Chromeleon 6.80 Service Pack 1 and later support Diagnostics functions for all pumps of the UltiMate 3000 system, except for the semipreparative HPG-3200P pump. These tests allow you to check the performance of certain pump components as well as the overall performance of the instrument. The Diagnostics Tool kit for the UltiMate 3000 pumps (part no. 6035.3000) includes all materials required for performing the tests. For the biocompatible pumps, an additional Adapter and Blind Nut kit (part no. 6037.3012) is required.

- 1. On the **Control** menu, select **Diagnostics**. (The **Control** menu is visible only when a control panel is open.)
- 2. The **Diagnostics** dialog box lists the available tests for all devices that are installed in the current timebase. Select a test for the pump. For information about how to run the tests, see the Chromeleon Help.

| To check the | Run the |
|---|----------------------|
| Pump system (especially the fitting connections) for leakage | General Leak Test |
| Pump valves and seals for leakage. The test helps you to identify the source for the leak. | Detailed Leak Test |
| Overall performance of the pump, that is the pressure pulsation and compressibility compensation | Performance Test |
| Backpressure caused by the filter frits in the inline filter of the outlet unit | Outlet Frit Test |
| Degasser vacuum. You can run this test for both the internal degasser of an LPG-3400 pump and the degasser of an SRD-3x00 Solvent Rack. | Degasser Vacuum Test |

If a test fails, check the Chromeleon Diagnostics Messages section for a short description of possible causes along with recommended courses of action (\rightarrow page 101).

5.6.3 Setting a Gradient Curve

For ramp gradients and multi-step gradients, you can specify linear or non-linear (curved) gradient profiles. The gradient profile (curve) is set in the program.

- 1. Create a program with the Program Wizard (\rightarrow page 65).
- 2. On the **Pump Options** page, select **Ramp** or **Multi-Step Gradient** on the **Gradient Type** list.
- 3. On the Flow Gradient Pump Options page, enter the desired curve (1 through 9) in the Curve field.

Curve 5, which is the default setting, is linear. Changes in composition of the delivered solvent over time are constant. Curves 1 through 4 are convex upward. Curves 6 through 9 are concave upward. For more information about gradient curves, see the Chromeleon Help.

5.6.4 Using the Digital Inputs and Outputs (Digital I/O)

You can specify in Chromeleon whether the pump controls the relays or whether the relays can be used in Chromeleon.

- 1. Open the **Commands** dialog box for the pump.
- For the relays R2 (LeftCamSyncOut), R3 (OperableOut), and R4 (RightCamSyncOut), the associated Enabled properties are available: Relay2Enabled, Relay3Enabled, and Relay4Enabled. To use a relay in Chromeleon, set the Enabled property for the relay to Yes. Set the Enabled property to No to have the pump control the relay for the associated signal.

If the relays are controlled by the pump, relays R2 and R4 provide the synchronization signal for the autosampler. The gradient running on the pump is synchronized with the **Inject** command of the autosampler.

Relay R2 can be controlled only by a DGP-3600 pump. With all other pumps, the relay can be used in Chromeleon in any case.

The **Operable Out** relay output (relay 3) closes when the pump is not ready to operate, that is, in case of an error or if the instrument is turned off.

- 3. On the **Relays** page of the pump properties, verify that the relays are enabled $(\rightarrow \text{ page 43})$. If the relay is not enabled, it will not be available in Chromeleon.
- **Tip:** The digital inputs are used, for example when a manual injection valve is connected to the pump (\rightarrow page 184).

For information about the pin assignment and the signal levels, see page 186.

5.6.5 Operational Qualification und Performance Qualification

Operational Qualification and Performance Qualification allow you to check and document the performance of the HPLC system. All materials required for performing qualification and detailed instructions are available on request.

5.6.6 Enabling the Double Flow Mode

Relevant only for HPG-3x00 pumps

In Double flow mode, both pump blocks are used for the delivery. In this way, the maximum allowed flow rate is doubled (\rightarrow page 156).

- 1. Start the **Server Configuration** program (\rightarrow page 36).
- 2. Right-click the pump in the timebase and select **Properties** on the menu.
- 3. On the **Head Type & Limits** page, select the **Double flow** check box (\rightarrow page 43).

5.7 Shutting Down the Pump

Observe the following precautions before interrupting the operation or before shipping the pump:

- Rinse out any solvents if necessary. If a buffer is used as a part of the mobile phase, flush the system with several volumes of methanol/water (50:50) before it is shut down. This will prevent salt buildup inside the unit.
- If operation is interrupted for more than 5 days, Thermo Fisher Scientific recommends removing the tubing from the peristaltic pump (→ Fig. 19, page 49). To remove the tubing, slightly press the lever to the right, remove the tubing, and release the lever. This will prevent that the tubing remains compressed and does not relax, thus blocking the wash solution.
- If operation is interrupted for more than one week, fill the pump with methanol or a similar alcohol, such as 2-propanol or ethanol. If the solvents in the pump are not miscible with water, use an appropriate intermediate solvent.
- Do not forget to empty the liquid reservoir of the rear seal wash system before shipping the pump (\rightarrow page 49).
- Rinse out buffers or solvents that form peroxide.
- Turn off the lamps in any UV or RF detectors that are connected to the pump. This will prevent evaporation of the solvents in the flow cell.
- Ship the unit only in the original shipping container and observe the packing instructions. Shipping the unit in any other packaging automatically voids the warranty. For more information, see the warranty statement in the terms of sale.

If you are running Chromeleon, you can set the pump and HPLC system into the standby mode or automate system shutdown.

Standby Program

A standby program sets the HPLC system into standby mode. The application can be reactivated very quickly afterward. The main program steps:

- The pump flow is automatically reduced at the end of the program.
- The temperature of all temperature-controlled modules in the system is reduced.

Shutdown Program

A shutdown program automates shutdown of the HPLC system. The main program steps are:

- The pump flow is automatically stopped at the end of the program.
- Certain system components and functions are turned off (for example, detector lamps, temperature control).

To create a standby or shutdown program

Select one of the following alternatives:

- Select and perform the operating commands and parameters from the **Commands** dialog box (→ page 62).
- Create and run a corresponding program to automate the process (\rightarrow page 65).
- Use the SmartShutdown Wizard to create and run the program (see below).

To create the program with the SmartShutdown wizard

- 1. To open the wizard, click **SmartShutdown** on the **Batch** menu.
- 2. Follow the instructions as they appear on each page of the wizard. For additional information about a page, click **Help**.
- 3. After you finish the wizard, Chromeleon
 - creates the program and saves it in the timebase for which you create the program.
 - opens the **Start Batch on** dialog box.

Select the program and click **Start** to run the program.

For more information about the SmartShutdown wizard, see the Chromeleon Help.

5.8 Routine and Preventive Maintenance Intervals

The pump is made of high-quality components and materials to minimize maintenance requirements. All surfaces are resistant to weak acids, alkali, and organic solvents. Nevertheless, immediately wipe up all liquids spilled onto the pump surface, using lint-free cloth or paper. If surfaces are exposed for longer periods, these liquids can cause damage.

Perform the maintenance procedures listed in the following table at regular intervals to ensure optimum performance and maximum uptime of the pump. The exact maintenance schedule for the pump will depend on a number of factors.

| Frequency | What you should do |
|-----------|--|
| Daily | Before you start operating the pump, inspect the fluid lines for air bubbles and degas the solvent. |
| | Check the fluid lines for indications of leakage. |
| | Check the fluid connections for indications of salt deposits. |
| | When using buffer solutions, flush the system thoroughly after use. Use a solvent that does not contain buffers or salts. |
| Regularly | Inspect the tubing for possible damage, such as cracks, nicks, cuts, or blockage. |
| | Check the liquid level in the liquid reservoir of the seal wash system. When using the internal reservoir, make sure that the level is always between the min. and max. markers on the label. |
| | Replace the liquid in the liquid reservoir of the seal wash system at regular intervals. <i>Internal liquid reservoir:</i> This is especially important when you use solvents with a high salt content. To avoid salt concentration in the reservoir and thus, to prevent damage to the piston seals, replace the liquid in the reservoir at least once a week. |
| | Inspect the filter frit in the inline filter for contaminants (\rightarrow page 112). Replace the filter frit if necessary (\rightarrow page 112). |
| | When using buffer solutions, Thermo Fisher Scientific recommends inspecting the pump for leakage at least once a month (\rightarrow page 117). |
| | Check the filter frits in the solvent lines for permeability. Replace the filter frits in regular intervals when using aqueous solvents. Aqueous solvents may contaminate the filters with algae and other microorganisms that deposit on the filter frits. Therefore, also replace the solvents at regular intervals. Rinse the reservoirs thoroughly before refilling them. |

| Frequency | What you should do | |
|-----------|--|-----------|
| Regularly | Check the drain tube connected to the drain port on the bottom right of the pump (\rightarrow Fig. 23, page 53). Verify that the tubing is unclogged and is routed below the drain port. Check the volume of the liquid in the waste container and empty as needed. | |
| Annually | Thermo Fisher Scientific recommends performing preventive maintenance on each pump annually. Preventive maintenance kits include all the required components: | |
| | Maintenance Kit for | Part No. |
| | ISO-3100A, LPG-3400A, LPG-3400M, DGP-3600A, DGP-3600M, HPG-3x00A, and HPG-3x00M | 6035.1961 |
| | LPG-3400AB, LPG-3400MB, DGP-3600AB, and DGP-3600MB | 6035.1963 |
| | HPG-3200P | 6035.1962 |
| | For information about the kits contents, see secti $(\rightarrow page 169)$. | on 11.3 |

Tip: Chromeleon supports functions for estimating the lifetime of consumables and diagnostic tests to check the performance of certain pump components $(\rightarrow \text{ pages 86 and 87}).$

6 Troubleshooting

6.1 Overview

The following features help you to identify and eliminate the source for problems that may occur during the operation of the pump.

Status Indicators

The status indicators on the front panel provide a quick visual check of the operational status of the pump. They indicate whether the pump is turned on, connected in Chromeleon, and operating properly (\rightarrow page 17).

Messages

If a fault or error is detected during the operation of the pump, a message is generated on the user interface. Check the Messages on the Front Panel Display section for recommended courses of action (\rightarrow page 96). If the pump is operated by Chromeleon, the message also appears in the Chromeleon Audit Trail.

Tip: For information about operating problems that might occur during the operation of an UltiMate 3000 system, see Operating Problems (\rightarrow page 103).

Diagnostics Tests

If the pump is connected in Chromeleon, Chromeleon provides several diagnostic tests allowing you to check the performance of certain pump components (\rightarrow page 87). If a test fails, check the Chromeleon Diagnostics Messages section for a short description of possible causes along with recommended courses of action (\rightarrow page 101).

Tip: To test the pump for leakage when the pump is not connected in Chromeleon, see section 7.8 (\rightarrow page 133).

If you are unable to eliminate a problem following the instructions given here, contact Thermo Fisher Scientific Service for Dionex HPLC Products.

6.2 Messages on the Front Panel Display

Each time a fault or error occurs during the operation of the pump, the **Status** LED on the front panel changes to red and a message appears on the front panel display. In this case, the **Prev**, **Next**, and **Clear** keys appear on the navigation bar.

| То | Select |
|---------------------------------------|--------|
| Return to the previous message. | Prev |
| Proceed to the next message. | Next |
| Remove all messages from the display. | Clear |

These keys are active also when the pump is connected in Chromeleon.

When the pump is operated with Chromeleon

- The error is also logged in the Chromeleon Audit Trail,. For some messages, the name specified on the **Devices** page in the Server Configuration program (→ page 39) may appear in front of the actual message text, for example, MicroPump.
- Messages on the front panel display can be removed also via the **ClearDisplayError** command in Chromeleon.

The following table lists pump-related messages and suggests appropriate remedial actions. In addition to the messages in the table, other messages may appear. In this case, note the exact wording of the message and contact Thermo Fisher Scientific Service for Dionex HPLC Products if you are unable to eliminate the problem.

| Message | Remedial Action |
|---|--|
| Abnormal drive current for xx seconds since (= time). | Inspect the capillary from the working head to the equilibration head for indications of blockage. Replace the capillary if necessary. Check the valve for permeability. |
| Camshaft index too early or too late. | Turn the pump off and on again by pressing the power switch on the rear of the pump. |
| Camshaft sensor always alight. | Turn the pump off and on again by pressing the power switch on the rear of the pump. |
| Camshaft sensor missing or dark. | Turn the pump off and on again by pressing the power switch on the rear of the pump |
| Degasser malfunction. | <i>LPG-3400</i> The vacuum level monitoring function of the degasser in an LPG-3400 pump recognized insufficient vacuum. Turn the pump off and on again by pressing the power switch on the rear of the pump. Check the degasser vacuum in Chromeleon (Commands dialog box). After about 1 minute, the setting should change from NotOk to OK. Inspect the degassing module for indications of leakage if the degasser does not reach the appropriate operating vacuum. |

| Message | Remedial Action |
|---|---|
| Degasser malfunction. (Cont'd) | All pumps except LPG-3400 The vacuum level monitoring function of the degasser in the SRD- 3x00 Solvent Rack recognized insufficient vacuum. If an SRD-3x00 Solvent Rack with integrated degasser is connected to the pump, turn the rack off and on again via the standby button. |
| Drive current too large. Camshaft xx°. | <i>Note: When this message appears, the pump continues running.</i> Contact Service. |
| Excessive drive current. Camshaft xx°. | Note: When this message appears, the pump continues running. The flow path before the system pressure sensor is blocked. Make sure that neither the capillary from the working cylinder to the equilibration cylinder nor the capillary to the outlet unit is blocked. Replace the capillaries if necessary. Verify that the valve cartridge in the check valve is installed in the direction of flow (\rightarrow page 112). Inspect the purge valve for indications of blockage (\rightarrow Fig. 9, page 23). |
| Invalid flow value. | The value you entered is not valid for this pump version configuration. Enter a flow value that it is within the permissible range. |
| Invalid parameter. | Enter a valid parameter. |
| Invalid partial flow setting. | The entered partial flow setting is not valid. Enter a partial flow setting that is within the permissible range and is logically correct. |
| Leak detected. | The leak sensor has reported a leak. Find and eliminate the source for the leak. Dry the leak sensor and tray (\rightarrow page 110). |
| Leak detected. Flow stopped. | The leak sensor has reported a leak. The flow is automatically turned off. Find and eliminate the source for the leak. Dry the leak sensor and tray (\rightarrow page 110). |
| Left hand pump block carryover pressure is too high. | Make sure that neither the capillary from the working cylinder to the equilibration cylinder nor the capillary to the outlet unit is blocked. Replace the capillaries if necessary. Inspect the purge valve (→ Fig. 9, page 23) for indications of blockage. |
| Motor malfunction. | Turn the pump off and on again by pressing the power switch on the rear of the pump. |
| Motor position error. The motor is overloaded. | Inspect the system for indications of blockage. This message can also appear if the upper pressure limit is exceeded (→ page 100). |
| Parameter could not be changed while pump is running. | You tried to change a parameter in a program while the pump or a program was running. Stop the pump or program and try again. |

| Message | Remedial Action |
|---|---|
| Possible piston seal leak detected. | More drops than specified under RearSealLeakage > Limit in Chromeleon reach the liquid reservoir. Visually inspect the piston seals for leakage (\rightarrow page 117). Replace the piston seals and support rings as necessary (\rightarrow page 119) or increase the default value for the piston leak in Chromeleon (\rightarrow RearSealLeakage > Limit , page 82). |
| Pressure fallen below lower limit. | The solvent supply is used up. Refill the solvent; purge the system $(\rightarrow page 54)$. There are air bubbles in solvent line. Check the filter frits; purge the system (\rightarrow page 54). The solvent emits gas when mixing. Degas the solvent and check the degasser. There is a leak in the system. Find and eliminate the source for the leak; tighten loose fitting connections. One of the check valves is defective. Clean or replace the valve (\rightarrow page 112). Purge the system as necessary (\rightarrow page 54). The lower pressure limit cannot be reached for the specified flow. The main task of the lower pressure limit is to monitor the system for leakage. Leakage may occur at the fittings and or screw joints, the high-pressure switching valve, or the piston seals. |
| Rear seal wash system has run out of wash solution. | Rear seal washing is enabled and the peristaltic pump is running, but no liquid reaches the seal-washing detector. Make sure the liquid reservoir contains sufficient liquid. Verify that the peristaltic tubing is connected correctly (→ Fig. 7, page 22). Verify that the peristaltic tubing is not blocked or drawing air. Replace the tubing as necessary. Clean the detector on the liquid reservoir. Make sure that the seal washing liquid is sufficiently conductive (→ page 49). |
| Relay 2 is configured for inject synchronization. Please change pump configuration. | You tried to use relay 2 as a standard relay in a program, although it has been set to synchronize the gradient with the Inject of the autosampler. In Chromeleon, set the EnableLeftCamSyncOut property to Off or use a different relay. |
| Relay 3 is configured as 'Operable' Out. Please change pump configuration. | You tried to use relay 3 as a standard relay in a program, although it has been configured as Operable Out. In Chromeleon, set the EnableOperableOut property to Off or use different relay. |
| Relay 4 is configured for inject synchronization. Please change pump configuration. | You tried to use relay 4 as a standard relay in a program, although it has been set to synchronize the gradient with the Inject of the autosampler. In Chromeleon, set the EnableRightCamSyncOut property to Off or use a different relay. |
| Message | Remedial Action |
|--|---|
| Right hand pump block carryover pressure is too high. | Make sure that neither the capillary from the working cylinder to the equilibration cylinder nor the capillary to the outlet unit is blocked. Replace the capillaries if necessary. Inspect the purge valve for indications of blockage (\rightarrow Fig. 9, page 23). |
| Solvent rack leak detected. | The leak sensor in the Solvent Rack has reported a leak. There is a leak in the system or a fluid connection is loose. Find and eliminate the source for the leak. Tighten leaking connection and dry the leak sensor (\rightarrow Solvent Rack Manual). |
| The maximum purge pressure was exceeded. | After you have initiated purging in Chromeleon (Purge property set to On), a pressure of more than 50 bar built up. |
| <i>Or</i> The pressure during purge exceeded 50 bar. | Verify that the purge valve is open (\rightarrow Fig. 9, page 23). Loosen the purge valve screw if necessary. If the message appears again, check the flow path for indications of blockage (see "The system pressure exceeded the safety limit"). |
| The pressure sensors have detected a disturbance. | The error can be due to large amounts of air in the pump heads, for example, during initial installation. |
| | Purge the pump blocks (\rightarrow page 54). |
| The pressure in the left (right) hand working cylinder | Inspect the system for indications of blockage (see "The system pressure exceeded the safety limit"). |
| exceeded the safety limit. | Especially inspect the capillary from the working head to the equilibration head. |
| | Verify that the valve cartridge in the check valve (outlet) is installed in the direction of flow (\rightarrow page 112). |
| The rear seal leak count is xx (counted drops) and has | More drops than specified under RearSealLeakage > Limit in Chromeleon reach the liquid reservoir. |
| exceeded the limit of yy (leak detection threshold) | Visually inspect the piston seals for leakage (\rightarrow page 117). |
| detection uneshold). | Replace the piston seals and support rings as necessary $(\rightarrow \text{ page 119})$ or increase the default value for the rear seal leak limit in Chromeleon ($\rightarrow \text{RearSealLeakage} > \text{Limit}$, page 82). |
| The rear seal leak sensor is malfunctioning. | The detector on the liquid reservoir is very dirty. Clean the electrodes of the detector, using water or solvent. If the message appears again, contact Service. |
| The system pressure exceeded the safety limit. | Open the purge screw to check if the flow path behind the outlet unit is blocked (\rightarrow Fig. 9, page 23). |
| | If it is - The column may be contaminated. Rinse or replace the column. - The sampler may be blocked. Eliminate the source for the blockage. If it is not Inspect the inline filter in the outlet unit for indications of |
| | blockage. - Replace the filter frit (\rightarrow page 111). |
| This function cannot be adjusted by the user. | Only authorized service personnel is allowed to change this parameter. |

| Message | Remedial Action |
|--------------------------------|---|
| Upper pressure limit exceeded. | See "The system pressure exceeded the safety limit." If the problem occurs due to column ageing, it may be sufficient to increase the setting for the upper pressure limit. |

When the pump is operated with Chromeleon and if communication between Chromeleon and the pump cannot be established, related messages may appear in the Chromeleon Audit Trail.

| Message | Remedial Action | |
|--|--|--|
| LPG-3400@USB-01610103 - Device not found on the USB. | The USB connection between the pump and the Chromeleon server may be interrupted. Check the USB connection. | |
| | The power supply to the pump may be interrupted. Check the power supply connection of the pump. | |
| Error opening LPG-3400 @USB-01610103 – The System | The USB connection between the pump and the Chromeleon server may be interrupted. Check the USB connection. | |
| cannot find the file specified | The power supply to the pump may be interrupted. Check the power supply connection of the pump. | |
| Error issuing control request to LPG-3400@USB-01610103 | The USB connection between the pump and the Chromeleon server may be interrupted. Check the USB connection. | |
| | The power supply to the pump may be interrupted. Check the power supply connection of the pump. | |
| | Remove the pump specified in the message from the server configuration or else, select a different pump from the list of available pumps in the server configuration program (via Properties and Browse). | |
| Error reading from LPG-3400 @USB-01610103 Data error (cyclic redundancy check) | Check the USB connection. The connection to the next hub must not exceed 5 m. The overall connection length, including the hub connections must not exceed 30 m. | |
| | Replace any defective USB cable or hub. | |
| Error reading from LPG-3400 @USB-01610103 | The USB connection between the pump and the Chromeleon server may be interrupted. Check the USB connection. | |
| | The power supply to the pump may be interrupted. Check the power supply connection of the pump. | |

6.3 Chromeleon Diagnostics Messages

If the pump fails a diagnostics test, perform the instructions given here and repeat the test. If the pump still fails the test, contact Thermo Fisher Scientific Service for Dionex HPLC Products.

Degasser Vacuum Test failed

| Possible Cause | Remedial Action |
|---|---|
| The degasser module did not reach the appropriate operating vacuum. | Check the tubing of the vacuum system and the degasser module. Repeat the test. If the test fails again, the degasser module may be defective. Contact Service. |

Detailed Leak Test failed

| Possible Cause | Remedial Action | |
|-------------------------------|--|--|
| The check valve has a leak. | Replace the outlet check valve (\rightarrow page 112) and repeat the test. | |
| | Replace the inlet check valve (\rightarrow page 112) and repeat the test. | |
| The piston seals are leaking. | Replace the piston seals in the working and equilibration heads $(\rightarrow page 119)$. | |

General Leak Test failed

| Possible Cause | Remedial Action |
|---|---|
| The pressure drop has been too high. | Check the blind put on the nump outlet for tightness |
| The pressure level has not been reached in time. | Visually inspect all fluidic connections for liquid droplets. |
| The initial pressure has not been reached within the first minute. | Consider replacing the inlet check valve (\rightarrow page 112). |
| The pressure drop during relaxation has been too high. The pump has a leak. | If you suit suspect there is a leak, full the detailed leak test. |

Outlet Frit Test failed

| Possible Cause | Remedial Action |
|--|--|
| The backpressure has exceeded the limit. The filter frit in the inline filter of the outlet unit may be blocked. | Replace the filter frit (\rightarrow page 112). |

Performance Test failed

| Test Result | Possible Cause | Remedial Action |
|---|---|---|
| The pressure/% compression was outside specification! | Eluent may not have been degassed correctly. | Degas the eluent. |
| | The wrong eluent may have been chosen in test step 1 (Prepare Pump). | Repeat the test. Make sure that the correct eluent is chosen. |
| | There may be air bubbles in the pump head. | Purge the pump (\rightarrow page 54). |
| | The check valves may not be working correctly. | Consider replacing the check values $(\rightarrow \text{ page 112}).$ |
| | Seals may not be tight. | Consider replacing the piston seals $(\rightarrow page 119).$ |
| | Backpressure may have fluctuated due to problems with the restrictor. The restrictor may be blocked. | Consider replacing the restrictor. |
| The pressure was outside specification. | The pressure was outside the upper and lower pressure limits (that is, $p < 50$ bar and $p > 400$ bar). | Repeat the test. Make sure that the pressure setting is correct. |
| The % compression value was over the limit! | The wrong eluent may have been chosen in test step 1 (Prepare Pump). | Repeat the test. Make sure that the correct eluent is chosen. |
| | The flow may have been too high. | Repeat the test with reduced flow. |
| The % compression value was below the limit! | It is likely that the wrong eluent was chosen in test step 1 (Prepare Pump). | Repeat the test. Make sure that the correct eluent is chosen. |
| | The flow may have been too low. | Repeat the test with increased flow. |
| Pressure ripple was too high. | There may be a mechanical fault. | Repeat the test. If the test fails again, contact Service. |

6.4 Operating Problems

The following table provides information about common operating problems that might occur with an UltiMate 3000 system and lists probable causes, as well as remedial actions. For more information, also see the manuals for the other modules of the UltiMate 3000 system.



If the pump is part of an UltiMate 3000 Proteomics MDLC system for nano or capillary HPLC, also see the information in the UltiMate 3000 Proteomics MDLC system manual.

| Problem | Probable Cause | Remedial Action |
|--|--|---|
| No information appears on the front panel display. | The instrument is not connected to the mains. | Connect the power cord. |
| | The power is turned off. | Turn on the power. |
| | The instrument is in standby mode. | Press the Standby key on the front panel. |
| | The screen brightness or contrast is not adjusted correctly. | Adjust the brightness or contrast on the Configuration menu or in Chromeleon. |
| | The fuses blow. | Replace the fuses $(\rightarrow page 131).$ |
| | Replacement fuse blows. | Contact Service. |
| | An error occurred in the electronic system. | Contact Service. |
| The instrument does not work correctly when controlled by Chromeleon. | There is no connection between the instrument and the Chromeleon computer. | Check the USB cable and connection to the computer. |
| | The USB port on the computer is not ready for operation. | Check the USB port on the computer. |
| The system has very high backpressure. | One or more capillaries in the system are blocked or damaged by bending. | Check the capillaries in the system, remove the blockage, or replace the capillaries. |
| The system has very high backpressure at the column and pump. | The column is contaminated or blocked. | Rinse or replace the column. |
| The system has very high backpressure at the pump (the pressure at the column is normal). | The filter frit on the high-pressure side is dirty. | Replace the filter frit (→ page 111). |
| | The filter frit installed on the outlet block is not the appropriate choice for the pump type. | Make sure that the appropriate filter frit is installed (\rightarrow page 23). Replace the filter frit if necessary (\rightarrow page 111). |

| Problem | Probable Cause | Remedial Action |
|--|---|--|
| High baseline drift | The column is contaminated. | Clean or replace the column. |
| | The system is not sufficiently equilibrated. | Flush the system until equilibration. |
| | The solvent is degraded. | Use fresh solvent and check the eluent filter frits. In aqueous solvents, growth of microorganisms is possible. |
| | The environmental conditions are unstable. | Make sure that the temperature and the humidity are constant. Avoid draft. |
| | The detector has not yet reached the operating temperature. | Allow the full detector warm- up time. |
| | The solvent is delivered in circles. | Direct the solvent to waste. |
| | The flow cell is dirty. | Clean the flow cell. Replace the flow cell if necessary $(\rightarrow Detector Manual)$. |
| | The lamp is too old. | Replace the detector lamp $(\rightarrow Detector Manual).$ |
| High noise level, pulsation | There are pressure fluctuations from the pump. | Purge the pump (\rightarrow page 54). |
| | There are air bubbles in the system. | Purge the pump (\rightarrow page 54). |
| | The solvent is degraded. | Use fresh solvent. |
| | The detector is defective. | Contact Service. |
| | The lamp is too old. | Replace the detector lamp $(\rightarrow Detector Manual).$ |
| | The wavelength is wrong. | Select an appropriate wavelength. |
| | The Step is too small (for example, 0.1 sec.). | Increase the Step (for example, to 0.5 to 1 sec.). |
| Pressure pulsation or inconstant pressure | The solvent is not degassed sufficiently. | Degas the solvent. |
| | The solvent is degraded. | Use fresh solvent. |
| | There are pressure fluctuations from the pump. | Purge the pump (\rightarrow page 54). |
| | There are air bubbles in the system. | Purge the pump (\rightarrow page 54). |

| Problem | Probable Cause | Remedial Action |
|---|--|---|
| Pressure pulsation or inconstant pressure (Cont'd) | The inlet valve or outlet valve on the working head is dirty. | Clean or replace the valve as necessary (\rightarrow page 112). |
| | The filter frit on the high- pressure side is dirty. | Replace the filter frit $(\rightarrow page 111).$ |
| Peak Broadening, increased dead time | The capillary to the detector is too long or the inner diameter of the capillary is too large. | Change the capillary. |
| | The filter frits on the solvent lines are clogged. | Check the filter for permeability. Replace the filter frit if necessary $(\rightarrow page 47)$. |
| | The capillaries are clogged. | Replace the capillaries. |
| | The sample loop is clogged. | Replace the sample loop $(\rightarrow Autosampler Manual).$ |
| | The proportioning valve is defective. | Contact Service. |
| | The column is overloaded or contaminated. | Clean or replace the column. |
| | The solvent is degraded. | Use fresh solvent. |
| Additional peaks appear in the injection peak. | With gradients, the equilibration time after the flush cycle is too short or the dead volume is too high. | Increase the equilibration time or eliminate possibly existing dead volume. |
| Poor reproducibility | The autosampler draws air from the vial. | There is not enough amount of sample in the vial, the needle height setting is incorrect (\rightarrow <i>Autosampler</i> <i>Manual</i>), or there are too many replicates. |
| | There are air bubbles in the syringe. | Flush the syringe $(\rightarrow Autosampler Manual).$ |
| | The draw speed for the sample is too high. | Reduce the draw speed. |
| | The gas content of the sample is too high or saturated. | Reduce the draw speed. Degas the sample if possible. |
| | The needle is clogged or the needle tip is deformed. | Replace the needle $(\rightarrow Autosampler Manual).$ |
| | There are air bubbles in the flow path. | Perform a wash cycle $(\rightarrow Autosampler Manual)$. Non-degassed wash liquid is used. Degas the wash liquid $(\rightarrow Autosampler Manual)$. |

| Problem | Probable Cause | Remedial Action |
|-------------------------------|--|---|
| Poor reproducibility (Cont'd) | The autosampler, the injection valve, or the syringe valve is not tight. | Inspect the autosampler and the connections on the injection valve or syringe valve. Tighten leaking connections. It might be necessary to replace the injection valve or syringe valve. In this case, contact Service. |
| | Carry-over occurs in the system. | Flush the needle using an appropriate solvent. |
| | The capillary connections are not installed properly or they are not | Check and tighten the capillary connections. |
| | tight. | Exchange the needle port if necessary (\rightarrow <i>Autosampler Manual</i>). |
| | | Exchange the needle if necessary (\rightarrow Autosampler Manual). |
| | There are dead volumes in the capillary connections. | Replace the fittings. Make sure that the capillaries are connected correctly. |
| | The piston seals are not tight. | Check the value for rear seal leak counter in Chromeleon (RearSealLeakage > Limit). Replace the seals (\rightarrow page 119). |
| | There is air in the working head. | Purge the pump (\rightarrow page 54). |
| | There is pump pulsation. | Use degassed solvents. |
| | | Check the filter frits in the solvent lines for contamination and exchange as necessary. |
| | | The outlet check valve may be defective. Replace the valve (\rightarrow page 112). |
| | The gradient is irreproducible. | Change the gradient. Check the pump function and degassing. Check the filter frits in the solvent lines for contamination and exchange as necessary. |

| Problem | Probable Cause | Remedial Action |
|-------------------------------|--|---|
| Poor reproducibility (Cont'd) | The sample is unstable and decomposes. | Use new sample or change the conditions. |
| | The environmental conditions are unstable. | Make sure that the temperature and air humidity are constant. (Use a column thermostat.) Avoid draft. |
| | Contamination occurs somewhere in the system. | Flush the system using an appropriate solvent. |
| No flow | The system is leaking. | Find and eliminate the leak. |
| | The valve cartridges are not installed correctly in the check valve nuts (not in the direction of flow) or are defective. | Correctly install or replace the check valve cartridges (→ page 112). |
| | There is air in the solvent or in the pump heads. | Purge the pump (\rightarrow page 54) or check the degasser. |
| | The inline filter of the pump is clogged. | Replace the filter frit $(\rightarrow page 111).$ |
| Flow fluctuation | The inlet path is blocked. | Check the inlet lines, filter of the pump, proportioning valve etc. for signs of blockage. |
| | There is air in the inlet path. | Purge the (\rightarrow page 54) and check the degasser. |
| | The check valves are dirty or defective. | Clean or replace the check valves (\rightarrow page 112). |
| | The piston does not contact the magnet holder. | Check whether the piston is installed correctly. Reinstall the piston (\rightarrow page 125). |
| | The piston seals are not tight. | Check the value for rear seal leak counter in Chromeleon (RearSealLeakage > Limit). Replace the seals (→ page 119). |
| Poor degassing | There is a leak in the capillaries or solvent lines or there are loose connections. | Inspect the capillary and solvent tube connections for leakage; tighten loose fitting connections. |
| | The degasser is not working properly. | Inspect the degassing chamber for indications for leakage. Check the degasser vacuum. |
| | The flow rate is too high. | Reduce the flow rate. |

| Problem | Probable Cause | Remedial Action |
|---|---|--|
| Reproducible ghost peaks in the chromatogram. | The degassing channels are contaminated. | Rinse the degassing channels $(\rightarrow \text{ page } 132).$ |
| | The solvents are degraded or dirty or their purity is insufficient. | Use fresh and appropriate solvents. |
| | Contamination occurs somewhere in the system. | Flush the system using an appropriate solvent. |
| The rear seal wash system is leaking. | The silicone tubing is bent or blocked. | Verify that the silicone tubing is installed properly $(\rightarrow page 22).$ |

7 Service

7.1 General Notes

The following sections describe all procedures for the pump that the user may perform. All other maintenance and service procedures must be performed by Thermo Fisher Scientific service personnel.

Observe the following precautions:

- Observe all warning notes when carrying out maintenance or repair work.
- Keep in mind that the fluidic components of the system may be filled with toxic solvents. Before starting maintenance or repair procedures, rinse toxic solvents from the instrument and put on protective clothing.
- Use original Dionex spare parts only. Substituting non-Dionex parts or using non-Dionex accessories may impair the performance of the instrument, thereby voiding the product warranty. For more information, see the warranty statement in the terms of sale.
- Before returning any instrument for repair, contact Thermo Fisher Scientific Service for Dionex HPLC Products. An RMA (Return Material Authorization) number is required to track your instrument. Always use the original packaging and observe the packing instructions when shipping the pump. Shipping the pump in anything other than the original packaging voids the warranty. For more information, see the warranty statement in the terms of sale.

If the original shipping container is not available, appropriate shipping containers and packing material can be ordered from the Thermo Fisher Scientific sales organization for Dionex HPLC Products. The packing instructions are included in the "Installation and Qualification Documents for Chromatography Instruments" binder and are available on request.

For instructions on shutting down the pump, see page 90.

Tip: Do not forget to test the pump for leakage after you have carried out maintenance or repair work on the fluid connections (\rightarrow page 133).

7.2 Drying the Leak Sensor

The leak sensor is installed inside the pump and reports a leak when liquid collects in the drip tray under the fluid connections. Find and eliminate the course for the leak and dry the sensor.



Fig. 39: Leak sensor

- 1. Wait until the pressure is down to zero.
- 2. Turn off the pump.
- 3. Inspect the pump heads for signs of leakage. Tighten or replace leaking connections.
- 4. Dry the components. If the pump heads continue to show signs of leakage, replace the piston seals (\rightarrow page 119).
- 5. With a cloth or tissue, absorb all liquid that has collected in the tray. Make sure that you do not bend or damage the sensor.
- 6. Allow the sensor to adjust to the ambient temperature for a few minutes.
- 7. Turn on the pump.
- 8. If no errors are reported after turning on the pump, operation can be resumed.
- 9. Do not forget to test the pump for leakage after you have carried out maintenance or repair work on the fluid connections (\rightarrow page 133).
- **I** Tip: If the sensor is not dry, the Status LED remains red. If a message appeared on the front panel display, select Clear on the navigation bar to remove the message from the display.

When the pump is operated with Chromeleon

If you eliminate the leak and dry the sensor *before* the **LeakDelay** has expired, the LED is green again. If you eliminate the leak and dry sensor *after* the **LeakDelay** has expired, the LED becomes green when you turn on again the pump flow (for a DGP-3600, both pump flows).

7.3 Filter Holder and Filter Frit in the Inline Filter

The outlet unit is fitted with an inline filter. The inline filter consists of the filter holder and the filter frit.



Inline Filter (= Filter holder with filter frit)



Filter holder

Filter frit

Fig. 40: Outlet unit with inline filter

| Description | Part No. |
|--|-----------|
| Filter holder for stainless steel pump | 6030.2103 |
| Filter holder for biocompatible pump | 6030.2112 |
| 10 filter frits, stainless steel, porosity: 0.5μm for the LPG-3400M and DGP-3600M pumps | 6000.0045 |
| 10 filter frits, titanium, porosity: 2 μm for LPG-3400MB and DGP-3600MB pumps | 6268.0036 |
| 10 filter frits, titanium, porostity: 10 μm for LPG-3400AB and DGP-3600AB pumps | 6268.0032 |
| 10 filter frits, stainless steel, porosity: 10µm for ISO-3100A, LPG-3400A, DGP-3600A, and HPG-3x00 pumps | 6268.0031 |

i Tips: The accessories kit for some pump types includes filter frits with different porosities. Therefore, make sure not to confuse the frit types.

When using filter frits with a porosity of 0.5 μ m, Thermo Fisher Scientific recommends installing PEEK filter frits for pH values < 4.

7.3.1 Checking the Permeability of the Filter Frit

Check the permeability of the filter frit at regular intervals.

- 1. Disconnect the capillary from the pump outlet.
- 2. Have the pump deliver water at a flow rate of 2 mL/min. When the outlet is open, the pressure should be less than 5 bar.
- 3. *Optional* Replace the filter frit as necessary (\rightarrow section 7.3.2).
- 4. Reconnect the capillary to the pump outlet.
- 5. Test the pump for leakage (\rightarrow page 133).
- **Tips:** In Chromeleon, you can check the permeability of the filter frit by running the **Outlet Frit Test** (\rightarrow page 87).

After you have replaced the filter frit, update the related service information in Chromeleon. To do so, perform the **OutletFritChanged** command.

7.3.2 Exchanging the Filter Holder or Filter Frit

- 1. Disconnect the capillary from the pump outlet.
- 2. Remove the filter holder from the outlet unit, using the open-end wrench from the pump accessories kit. The filter frit is in the filter holder.
- 3. Remove the frit from the holder.
- 4. Insert the new frit into the holder. Wet the frit on the side that goes into the holder, for example, with Isopropanol. This keeps the frit in the holder when you tighten the assembly onto the outlet unit.
- 5. Tighten the filter holder onto the outlet unit hand-tight.
- 6. Reconnect the capillary to the pump outlet.
- 7. Test the pump for leakage (\rightarrow page 133).
- **Tip:** After you have replaced the filter frit, update the related service information in Chromeleon. To do so, perform the **OutletFritChanged** command.

7.4 Replacing the Check Valves

The working head contains two double check valves: an inlet valve and an outlet valve. The inlet valve is mounted on the bottom side of pump head; the outlet valve is mounted on its topside. When removed, you can recognize the inlet valve nut by the inner thread that ends in a planar surface, while the outlet valve nut has a conical cavity for the capillary connection.

| Beschreibung | BestNr. |
|---|-------------------------------------|
| Valve cartridge (sapphire) | 6035.2300 |
| Check valves (complete): inlet and outlet check valve nuts for Analytical and micro pumps (A and M pump versions) Biocompatible pumps (AB and MB pump versions) Semipreparative pump (HPG-3200P) | 6035.1965 6037.1964 6035.1966 |

- 1. If necessary, rinse the pump to remove toxic solutions.
- 2. Set the pump flow rate to 0. Wait until the pressure in the system is down to zero.
- 3. Disconnect the tubing connections from the inlet and outlet check valves.
- 4. Loosen both check valve nuts, using a 13 mm or 1/2" wrench, and remove them from the pump head.
- 5. Hold the nuts over your hand and turn them upside down, allowing the cartridges into your hand.
- 6. Insert a new cartridge into the check valve nut. Make sure that you insert the cartridge in the direction of solvent flow, as indicated by the arrow on the cartridge. Screw the check valve into the pump head and tighten the check valve
- 7. Reconnect the tubing connections to the inlet and outlet check valve nuts. On the outlet check valve, tighten the tubing hand-tight, and then tighten it an additional one-quarter turn, using a wrench. When connecting the tubing to the inlet check valve, take care to avoid cross-threading. (If the valve leaks, you may tighten the fitting a little more. Do not overtighten; this will crush the cartridge.)
- 8. To prevent contaminants from entering the HPLC system, thoroughly rinse the pump (using at least 30 mL HPLC-grade water or purely organic solution). Loosen the purge valve screw to prevent the rinsing liquid from entering the HPLC system.
- 9. Test the pump for leakage (\rightarrow page 133). Tighten any leaking connections.
- **Tip**: After you have replaced the check valves, update the related service information in Chromeleon. To do so, perform the **CheckValveServiceDone** command.

7.5 Pistons and Piston Seals

Analytical, micro, and biocompatible pump versions



Fig. 41: Pump heads, pistons, and piston seals (here all pumps except HPG-3200P)

Depending on the pump version (analytical, micro, or biocompatible), the design of some parts may differ from the representation above (for part number information, see the following tables). However, the position of the parts is the same for all pump versions.

i Tip: HPG-3x00 pumps only

The working head is fitted with a pressure transducer.

| No. | Description | Part No. | | |
|-----|--|----------------------------------|------------------------------------|------------------------|
| | | ISO-3100A | LPG-3400A DGP-3600A | HPG-3x00A HPG-3x00M |
| 1 | Holder + piston support hemisphere | 6269.0040 | (piston support hemis | phere only) |
| 2 | Piston | | 6035.2240 | |
| 3 | Retainer | | Included in 6035.2020 ³ | * |
| 4 | Spring | | | |
| 5 | Rear seal wash body | | Included in 6035.2020 ³ | * |
| 6 | Ring seal (DR-8) | 2266.0082 (10 seals) | | |
| 7 | Support ring (piston seal) | 6025.2010A (Reversed Phase) | | |
| 8 | Piston seal (1/8") | <u>6025.2011A (Normal Phase)</u> | | |
| 9 | Equilibration head | | | |
| 10 | Valve cartridge (same as no. 13) | 6035.2300 | | |
| 11 | Inlet check valve nut | Included in 6035.1965 | | |
| 12 | Working head | | | |
| 13 | Valve cartridge (same as no. 10) | 6035.2300 | | |
| 14 | Outlet check valve nut | Included in 6035.1965 | | |
| 15 | Capillary from working head to equilibration head (U tube) | 6030.3015 | | |

| No. | Description | Part No. | | |
|-----|--|---|--------------------------|--------------------------|
| | | LPG-3400M DGP-3600M | LPG-3400MB DGP-3600MB | LPG-3400AB DGP-3600AB |
| 1 | Holder + piston support hemisphere | 6269.0040 (| piston support hemisp | ohere only) |
| 2 | Piston | | 6035.2240 | |
| 3 | Retainer | Inclue 6035.2 | ded in 2010** | Included in 6035.2020* |
| 4 | Spring | | | |
| 5 | Rear seal wash body | Included in Included in 6035.2010** 6035.2020* | | Included in 6035.2020* |
| 6 | Ring seal (DR-8) | 2266.0082 (10 seals) | | |
| 7 | Support ring (piston seal) | 6025.2010A (RP) | 6025.201 | 2 (RP) |
| 8 | Piston seal (1/8") | 6025.2011A (NP) | | |
| 9 | Equilibration head | | | |
| 10 | Valve cartridge (same as no. 13) | 6035.2300 | | |
| 11 | Inlet check valve nut | Included in 6035.1965 | Includ 6037. | ed in 1964 |
| 12 | Working head | | | |
| 13 | Valve cartridge (same as no. 10) | 6035.2300 | | |
| 14 | Outlet check valve nut | Included in 6035.1965 | Includ 6037. | ed in 1964 |
| 15 | Capillary from working head to equilibration head (U tube) | 6030.3015 | 6037.2 | 3015 |

Part no. 6035.2020 (rear seal wash system) contains 1 piston seal, rear seal wash body and retainer each.
 Part no. 6035.2010 (rear seal wash system) contains 1 piston seal, rear seal wash body and retainer each.

Semipreparative pump version



Fig. 42: Pump heads, pistons, and piston seals (here: HPG-3200P)

Semipreparative pump version

| No. | Description | Part No. |
|-----|---|-------------------------------------|
| 1 | Piston | 2267.0802A |
| 2 | Piston retainer | 6004.8104 |
| 3 | Piston bushing (5 bushings) | 2090.8083 |
| 4 | Retainer | Included in 6025.2210A* |
| 5 | Rear seal wash body | Included in 6025.2210A* |
| 6 | Ring seal (10 seals) | 2266.0082 |
| 7 | Support ring (piston seal) | 6030.9010 (für Reversed Phase) |
| 8 | Piston seal | $\int 6030.9011$ (für Normal Phase) |
| 9 | Capillary from working head to equilibration head (U tube) | 5030.3025 |
| 10 | Outlet check valve nut | Included in 6035.1966 |
| 11 | Valve cartridge (same as no. 14) | 6035.2300 |
| 12 | Working head with pressure transducer | |
| 13 | Inlet check valve nut | Included in 6035.1966 |
| 14 | Valve cartridge (same as no. 11) | 6035.2300 |
| 15 | Equilibration head (without pressure transducer) | |
| 16 | Connect the capillary from left equilibration head to Tee insert right equilibration head to Tee insert | Included in 6035.1969 |

* Part no. 6025.2210A (rear seal wash system) contains 1 piston seal, rear seal wash body and retainer each.

7.5.1 Visually Inspecting the Pump for Piston Seal Leakage

Each piston has two piston seals. The seals prevent solvent from leaking either into the rear seal wash system or over the piston and into the instrument. With the standard setup of the seal wash system (that is, the liquid reservoir is installed inside the pump), the detector on the liquid reservoir automatically checks the main piston seals for leakage (\rightarrow page 82). If rear seal washing is performed from an external liquid reservoir (\rightarrow page 83), the detector is not active.

In both cases, you can inspect the pump also visually for liquid leaks from the main piston seals.

Standard Setup with Internal Wash Liquid Reservoir

 Turn rear seal washing off and on again. Open the Commands dialog box for the pump in Chromeleon and set RearSealWashSystem first to Off and then to Automatic. —Or—

On the front panel display, select the **Preferences** menu and set **Rear seal wash** first to **Off** and then to **Automatic**.

The peristaltic pump starts pumping, delivers approximately 2 mL of liquid, and then stops.

2. Remove the silicone tubing from the pump head on the liquid reservoir (detector).



Fig. 43: Tubing connections on the internal liquid reservoir

- 3. Remove some of the liquid by shaking the tube.
- 4. Reconnect the tubing to the liquid reservoir.
- 5. Set the flow rate. Arrange the system in such a way that approximately 300 bar of backpressure is generated.

6. To evaluate possible leakage, observe the air/liquid level in the silicone tube.

I Tip: The peristaltic pump will start a new washing cycle after one hour. Therefore, be sure to finish your observation before that time.

7. If the level remains unchanged, the piston seals seal tightly. A rising or falling level indicates leakage from one or more of the main piston seals. In this case, replace all piston seals and the support rings (\rightarrow page 123).

Alternative Setup with External Wash Liquid Reservoir

Bad retention time precision may indicate piston seal leakage. Observe the open end of the waste line. If liquid exits the waste line during the intervals indicates liquid leaks from the main piston seals. In this case, replace all piston seals and the support rings (\rightarrow page 123).

7.5.2 Replacing the Piston Seals

A defective piston seal allows leakage past the piston. This may cause unstable flow rates and baseline noise; in addition, it may make it difficult to prime the pump.

The seal replacement procedure consists of:

- 1. Removing the pump head and piston (\rightarrow page 120)
- 2. Removing the piston seals and support ring (\rightarrow page 123)
- 3. Cleaning the pistons (\rightarrow page 124)
- 4. Installing the piston, piston seals, support ring, and pump head (\rightarrow page 125)
- 5. After you have replaced the piston seals, observe the recommendations on page 130.
- **I** Tips: As standard, all pumps are equipped with piston seals made of UHMW-PE (RP). Keep in mind that using chloroform, trichlorobenzene, methylene chloride, tetrahydrofuran, or toluene as solvents chemically damages the UHMW-PE seals. Chemical reactions may also occur when using tetrachloromethan, diethyl ether, ethyl ether, di-isopropyl ether, ketones, methylbenzene, methycyclohexane, and monochlorobenzene. If you use these solvents, contact the Thermo Fisher Scientific sales organization for Dionex HPLC Products.

As standard, the pumps are fitted with reversed phase piston seals. However, normal phase seals can be installed instead if required. Depending on the application, the abrasion may be slightly increased. Therefore, check the permeability of the filter frit at regular intervals (\rightarrow page 112).

For information about the installation procedure and the corresponding part numbers, see section 7.5.2.2 (\rightarrow page 123).

7.5.2.1 Removing the Pump Head and Piston

To reach the piston heads and pistons, it is not necessary to open the pump enclosure. Tilt the front cover upward. The pistons are freely accessible via the pump heads.

All pumps

- 1. If necessary, purge the pump to remove toxic solvents.
- 2. Set the pump flow rate to 0. Wait until the system pressure is down to zero.
- 3. Disconnect all fluid connections from the pump head.



Fig. 44: Installed pump heads

4. Loosen the Allen screws on the two pump head (\rightarrow Fig. 44) and carefully remove the pump head.



Fig. 45: Pump head removed

- 5. Remove the support ring. The support ring may still be in the removed pump head.
- 6. Continue with the steps that are appropriate for your pump.
 - For all pumps except HPG-3200P (\rightarrow page 121).
 - For HPG-3200P pumps (\rightarrow page 122).

All pumps except HPG-3200P

Carefully remove the rear seal wash body from the piston (→ Fig. 47), by pulling it toward you. To do so, use the extractor from the pump accessories kit (→ Fig. 46). Avoid tilting. This may cause damage to the piston.





Fig. 46: Extractor for rear seal wash body

Fig. 47: Extractor on rear seal wash body

8. Hold the piston as far as possible at its rear end. Move the piston, for example, upward, so that it is no longer in the center position.



Piston moved upward

Rear end of the piston

Fig. 48: Piston moved upward

- 9. Tilt the piston and remove it from the magnet holder (\rightarrow Fig. 41, no. 1, page 114).
- 10. Verify that the hemisphere (part no. 6269.0040), which supports the piston, remained in the holder. In rare cases, the hemisphere sticks to the piston and is removed with it. In this case, rub the hemisphere with a dry tissue. Apply some grease to the center of the holder (thin film). Insert the hemisphere with tweezers. The flat side must point toward you (\rightarrow Fig. 49).



HPG-3200P only

- 7. Carefully remove the rear seal wash body from the piston (\rightarrow Fig. 45), by pulling it toward you.
- 8. Remove the piston bushing from the piston, using the extractor from the pump accessories kit.



Fig. 50: Piston with piston bushing

Fig. 51: Removing the piston bushing

9. Loosen the piston retainer (\rightarrow Fig. 42, no. 2, page 116), using the slotted screwdriver from the pump accessories kit, and remove the piston from the enclosure.

7.5.2.2 Removing the Piston Seals and Support Ring

Each piston has two seals: a main piston seal (on the pump head) and a seal in the rear seal system.

- **I** Tip: Use a disassembled piston to remove the piston seal from the pump head. Do not use a sharp tool, such as tweezers. This will scratch the inside of the pump head; these scratches will prevent a proper seal and allow leakage when you install the new seal.
- 1. On the working head (with valve cartridges), use a blind nut (FS-8, part no. 6000.0044, for all pumps except biocompatible pumps; FS-8 biocompatible, part no. 6000.0144, for biocompatible pumps) to close the outlet valve.
- 2. On the equilibration head, close the holes with blind nuts.
- 3. Place a few drops of water on the main piston seal (for example, with a syringe). The seal is easier to remove when it is moist.
- 4. Push the piston into the pump head. The pressure loosens the seal. If you cannot remove the piston seal in this way, use an M4 screw (this is possible only for analytical pumps); for example, the one used for holding the pump heads. Insert the screw into the seal and remove the seal. This procedure destroys the piston seal.

M Important: When replacing the piston seal in the pump head, always replace the support ring as well. (For part number information, see the following table.) This is to prevent leakage.

Even if you do not want to replace the seal in the rear seal body, visually inspect the seal to verify that it is neither worn nor bent. To see the seal, you have to remove the retainer (see the next step below).

▲ Important: Lorsque vous remplacez le joint de piston de la tête de pompe, remplacez systématiquement la rondelle anti-extrusion. (Pour des informations sur les références des pièces, référez-vous au tableau cidessous.) Ceci est destiné à éviter toute fuite.

> Même si vous ne voulez pas remplacer le joint dans le corps d'arrièrejoint, inspectez visuellement le joint pour vérifier qu'il ni n'est porté ni est plié. Pour voir le joint, vous devez enlever l'arrêtoir (voir la prochaine étape ci-dessous).

| Description | Part No. |
|--|---|
| Piston seal/support ring kit (reversed phase) for - ISO-3100A, LPG-3400A and M, DGP-3600A and M, and HPG-3x00A and M - LPG-3400AB and MB, DGP-3600AB and MB - HPG-3200P | 6025.2010A* 6025.2012* 6030.9010* |
| Piston seal/support ring kit (normal phase) for - ISO-3100A, LPG-3400A and M, DGP-3600A and M, and HPG-3x00A and M - HPG-3200P | 6025.2011A* 6030.9011* |

* The kits include one support ring and two piston seals.

 To remove the piston seal that is installed in the rear seal body, remove the retainer on the back of the rear seal wash body first, using the extractor tool from the accessories kit (→ Fig. 46, page 121 and Fig. 52, page 125). Then, remove the seal. Use the piston to push the piston seal out of the rear seal body.

The seal does not only seal the rear seal wash system but also guides the piston. Therefore, always replace the seal when it is worn or bent.

7.5.2.3 Cleaning the Pistons

All pumps

- 1. Remove the pump head and the piston if necessary (\rightarrow page 120).
- 2. Carefully rinse the piston, and then rub it several times with a dry, lint-free tissue.
- 3. Inspect the piston for signs of damage. If the piston is scratched or scored, replace it.

HPG-3200P only

- 4. Apply some thin-bodied, resin-free oil (thin film) only to the metal part of the piston.
- **I** Tip: Use a lamp to find out whether the piston is dirty.

Sapphire piston

Hold the rear side of the piston into the light. Dirt particles will be enlarged by the refraction of the light.

Ceramic piston

Hold the piston into the light. While turning the piston, observe the reflection of the lamp.

7.5.2.4 Installing the Piston, Piston Seals, Support Ring, and Pump Head

- **Tip:** Thermo Fisher Scientific recommends that you always install new piston seals when you install a piston.
- 1. Assemble the rear seal wash system. If you did not disassemble the rear seal wash system, continue with the step 5 that is appropriate for your pump.
- 2. Insert the piston into the rear seal wash body (from the side opposite the retainer; \rightarrow Fig. 52).
- 3. Slide the new piston seal (mind the correct orientation) over the piston. (The piston is used only to facilitate centering during this process.) Insert the retainer into the rear seal body. Tighten the retainer finger-tight, and then use extractor to tighten the retainer gently until the seal snaps into place. Remove the piston so that the seal remains in the rear seal body.



Fig. 52: Assembling the rear seal wash system

- ▲ Important: When installing the piston seals, make sure that the spring side of the seals faces away from the pump enclosure (→ Fig. 41 and Fig. 42, no. 8).
- ▲ Important: Lorsque vous installez des joints de piston, assurez-vous que les côtés "ouverts" des joints, montrant le ressort spiral, soient dirigés vers l'avant de la tête de pompe (coté pression) (→ Fig. 41 et Fig. 42, no. 8).
- 4. Continue with the steps that are appropriate for your pump.
 - For all pumps except HPG-3200P (\rightarrow 126).
 - For HPG-3200P pumps (\rightarrow 128).

All pumps except HPG-3200P

5. Verify that the O-ring seal (part no. 6266.2202 for 5 seals) is installed on the rear seal wash body (→ Fig. 53). Slightly grease the O-ring seal and metal part of the rear seal wash body if necessary.



Fig. 53: Rear seal wash body with O-ring seal

- 6. Insert the piston into the rear seal wash body only that far that you can slide the support ring and piston seal onto the other piston end. Then, slide the support ring and ring seal onto the piston.
- 7. Verify that the uninstalled pump head has no seal installed (for example, the old seal) and that the white ring seal is on the rear seal wash body (\rightarrow Fig. 45, page 120).

Slightly push the rear seal wash body into the pump head until the piston seal is correctly in place. Avoid tilting and only push the rear seal wash body. To avoid damage to the piston, do not push from the piston end.

- **Tip**: To avoid damage to the piston seal during installation, Thermo Fisher Scientific recommends moistening the piston seal, piston, and pump head with HPLC-grade water directly before you install them.
- 8. Slide the spacing tool over the piston. Move the tool forward until it is stopped by the rear seal wash body. Remove the tool. The piston is now in the correct position for being installed in the pump head.



Fig. 54: Spacing tool

Use this side of the tool with all pumps *except* LPG-3400M(B) and DGP-3600m(B).

Use this side of the tool (deeper spacing) with the LPG-3400M(B) and DGP-3600M(B).

- 9. Verify that the magnet holder is free from contaminants and that the piston support hemisphere is in the holder. It is important that the piston does not contact the magnet holder immediately upon installation. Therefore, perform the Change right pump pistons or Change left pump pistons command on the Control menu (→ page 71) or perform the related command in Chromeleon to move the pump in the appropriate position. Then, slide the pump head together with the rear seal wash body and piston into the pump block and tighten the Allen screws in turn.
 - **Important:** If you do not insert the piston seal in the way described above, but into the disassembled pump head, it may get jammed and thus become useless.

The working head is installed on the left side of the pump block (inlet). You can identify the head by the two large holes for the inlet and outlet valves.

▲ Important: Si vous n'insérez pas le joint de piston de la manière décrite cidessus, mais dans la tête de pompe démontée, il peut se coincer de travers et ainsi devenir inutilisable.

> La tête de travail est installée sur le côté gauche du bloc de pompe (entrée). Vous pouvez reconnaître la tête aux deux grands taraudages destinés aux supports des clapets d'entrée et de sortie.

- 10. Have the pump deliver about 1 mL/min (analytical pumps) or $300 \,\mu$ L/min (micro pumps). The magnet holder moves forward until the piston and magnet holder come into contact (indicated by an audible click). If you replace pistons, two clicks must be heard. Only then, both pistons are in the correct positions.
- 11. Reinstall the capillary connections as described in the All Pumps section $(\rightarrow \text{ page 129})$.

HPG-3200P only

- 5. Install the piston in the pump block and tighten the piston retainer using the slotted screwdriver from the accessories kit.
- 6. Slide the piston bushing onto the piston. Be sure to install the bushing in the correct orientation (\rightarrow Fig. 55).

Slide the bushing onto the piston in this direction.



Fig. 55: Correct orientation of the piston bushing

- 7. Slide the complete rear seal wash body over the piston.
- 8. Slide the new support ring and the new piston seal onto the piston. Be sure to install the piston seal in the correct orientation. When the seal is installed correctly, the open side points toward the front.



Fig. 56: Rear seal wash system with seals and support ring

- Slide the ring seal (part no. 2266.0082; 10 seals) onto the rear seal wash body (→ Fig. 42, no. 6).
- 10. Verify that the uninstalled pump head has no seal installed (for example, the old seal).
- 11. Slide the pump head onto the piston (turn and push gently as necessary) until the second piston seal is correctly in place.

Important: If you do not insert the piston seal in the way described above, but into the disassembled pump head, it may get jammed and thus become useless.

The working head is installed on the left side of the pump block (inlet). You can identify the head by the two large boreholes for the inlet and outlet valves.

- Important: Si vous n'insérez pas le joint de piston de la manière décrite cidessus, mais dans la tête de pompe démontée, il peut se coincer de travers et ainsi devenir inutilisable.
 La tête de travail est installée sur le côté gauche du bloc de pompe (entrée). Vous pouvez reconnaître la tête aux deux grands taraudages destinés aux supports des clapets d'entrée et de sortie.
 Tip: To avoid damage to the piston seal during installation, Thermo Fisher Scientific recommends moistening the piston seal, piston, and pump head with HPLC-grade water directly before you install them.
- 12. Tighten the pump head with the Allen screws.
- 13. Reinstall the capillary connections as described in the All Pumps section $(\rightarrow page 129)$.

All pumps

14. Reinstall all capillary connections to the pump heads. When you install the solvent line on the inlet valve, take care to avoid cross-threading.

HPG-3x00M only

Due to limited space, it may be difficult to install the capillary from the left equilibration head to the outlet unit. Be careful not to bend the capillary on the equilibration head; this will prevent a proper seal and allow leakage. To facilitate the installation, follow the steps below:

- a) Move the fitting screws on the capillary as far back as possible.
- b) First, insert the capillary into the hole on the equilibration head. Then, slightly turn the capillary as necessary and insert the other end into the hole on the outlet unit. If you need to bend the capillary during the installation, be careful to bend the tubing only slightly and only at the position marked in Fig. 57. Only at these positions, the capillary is flexible enough to return to its original shape after being bent.



Fig. 57: Capillary from left equilibration head to outlet unit (HPG-3x00M)

c) Move the fitting screws back and tighten the capillary.

- 15. Attach the tubing of the rear seal wash system to the corresponding capillaries (\rightarrow Fig. 7, page 22).
- 16. Rinse the pump thoroughly to prevent contaminants from entering the HPLC system.
 - Rinse the pump with at least 30 mL of HPLC-grade water or pure organic solution.
 - Loosen the purge valve screw to prevent rinsing liquid from entering the HPLC system.
- 17. After replacing the piston seals, observe the recommendations on page 130.
- 18. Test the pump for leakage (\rightarrow page 133). Tighten any leaking connections.
- **Tip:** After you have exchanged the piston seals or pistons, update the related service information in Chromeleon. To do so, perform the **PistonsChanged** and **SealsChanged** commands.

7.5.2.5 Recommended Actions after Exchanging the Piston Seals

Observe the following recommendation when you have exchanged the piston seals.

- Always exchange the liquid in the reservoir of the rear seal wash system and rinse the fluidics.
- Allow new piston seals to run in.
 - a) Connect drain tubing to the purge outlet.
 - b) Open the purge valve.
 - c) Operate the pump for 15 minutes with isopropanol at a flow rate of 1 mL/min. Do not have the pump deliver in circles.
 - d) Close the purge valve.
 - e) On the pump outlet, install a flow resistance, for example, a capillary, that can generate approximately 300 bar at a flow rate of 1 mL/min.
 - f) Have the pump deliver isopropanol at 1 mL/min for another 30 minutes.
 - g) Inspect the filter frit in the outlet unit for permeability (\rightarrow page 112).
 - h) Uninstall the capillary and connect the pump to the system.

In rare cases, it may happen that new seals still show an increased leakage rate after several hours of operation. In this case, exchange the liquid in the reservoir of the rear seal wash system once again and rinse the fluidics.

- If leakage is observed with new piston seals, operate the pump for at least 2 hours at 350 bar to run in the seals. *Note*: In this case, tightening the pump head screws does not help.
- Never run the pump dry. Damage to the pistons or piston seals could result.

7.6 Replacing the Fuses

STOP

- **Warning:** Before replacing the fuses, turn off the pump. Be sure to disconnect the power cord from its source.
- Avertissement: Avant de remplacer les fusibles, arrêtez la pompe. Assurez-vous de bien débrancher le cordon d'alimentation de la source secteur.
- 1. Remove the fuse cartridge, using a small screwdriver (\rightarrow Fig. 58).



Fig. 58: Fuse cartridge

- 2. Replace the fuses with fuses of the appropriate rating. Thermo Fisher Scientific recommends always replacing both fuses.
- **Important:** Do not operate the pump with only one fuse. Use only the fuses indicated in the following table.
- **Important:** Ne faites pas fonctionner la pompe avec seulement un fusible. Utilisez uniquement les fusibles indiqués ci-dessous.

| Description | Part No. |
|--|---|
| Fuse, 2 A, slow-blow, 5 x 20 mm, 250V | Included in Fuses Kit, part no. 6030.9003 For information about which fuses are contained in the kit, see section 11.3 (\rightarrow page 169). |

- 3. Reinstall the fuse cartridge.
- 4. Reconnect the power cord to its source and turn on the pump.

7.7 Vacuum Degasser

To avoid contamination of the degasser

- Prepare fresh solvents at regular intervals.
- Rinse the degassing channels at regular intervals.
 - This is especially important for the channel that degases the aqueous solvent (for reversephase chromatography). Usually, it is sufficient to rinse all channels with organic solvent. Use a fresh bottle. When you use water and acetonitrile or methanol, it is usually sufficient to rinse the degassing channels once per week. However, adapt the rinsing intervals to the solvents in use. In persistent cases, for example reproducible ghost peaks in the chromatogram, follow the steps below:
 - 1. Install a backpressure capillary on the pump outlet. The capillary should be appropriate for generating a backpressure of 200 to 300 bar.
 - 2. Rinse the degassing channels for 1 hour with 20% nitrid acid at the flow rate normally used for your application. Use fresh HPLC-grade water.
 - 3. Rinse the degassing channels with fresh HPLC-grade water until the pH value is neutral.
 - 4. Rinse the degassing channels for 2 hours, using fresh HPLC-grade acetonitrile quality at the flow rate normally used for your application.
 - 5. Prepare fresh solvents for your application using new bottles, and connect them to the degassing channels.
 - 6. Install new filter frits on the solvent lines.
 - 7. Uninstall the backpressure capillary and reconnect the system as required by your application.
 - 8. Equilibrate the system.

In addition, observe the following

- Clean the solvent lines at regular intervals.
- Thermo Fisher Scientific advises against recycling the solvents. This may impair the degassing performance.
- When connecting the solvent lines, make sure that the connectors are free of contaminants. Even minute particles can allow air to enter the degasser, and thus reduce the unit's effectiveness.
- When replacing solvents, make sure that the solvents are miscible. Mix immiscible solvents with an intermediate solvent to replace them step-by-step.
- Thoroughly rinse the degasser with alcohol after operation. The alcohol does not need to be removed afterwards.
- For longer periods of inactivity and when using saliferous buffers (which may result in salt crystallization in the gas separation membrane, thereby impairing the degassing performance), rinse with de-ionized water followed by either methanol or ispropanol.

7.8 Testing the Pump for Leakage

After you have carried out any maintenance or repair work on the fluid system, test the pump for leakage.

I Tips: Perform the test for the two pumps of a DGP-3600 separately if necessary.

In Chromeleon, you can run **General Leak Test** (\rightarrow page 87) to test the pump for leakage.

1. Close the pump outlet with an appropriate blind nut.

| Description | Part No. |
|---|-----------|
| Blind nut (FS-8) for all pumps except biocompatible pumps | 6000.0044 |
| Blind nut (FS-8) for biocompatible pumps | 6000.0144 |

 In Chromeleon, set the upper pressure limit: <u>For pumps without flow splitter assignment</u>: Set Pressure > UpperLimit. <u>For pumps with flow splitter assignment</u>: Set MasterPressure > UpperLimit.

| Ритр Туре | Upper Pressure Limit | |
|----------------------------|----------------------|--|
| Analytical and micro pumps | 400 bar | |
| Semipreparative pumps | 100 bar | |

- 3. In Chromeleon, select a flow rate of, for example, 50 μ L/min.
- 4. Decrease the flow as soon as the pressure builds up (see table).

| Ритр Туре | Pressure typically |
|----------------------------|-------------------------|
| Analytical and micro pumps | between 100 and 200 bar |
| Semipreparative pumps | between 80 and 100 bar |

5. Have the pump deliver some μ L/min until the appropriate pressure has build up.

| Ритр Туре | Pressure Build Up |
|----------------------------|-------------------------|
| Analytical and micro pumps | between 100 and 200 bar |
| Semipreparative pumps | between 80 and 100 bar |

6. When this pressure has been built up and when the pump delivers a master flow of $1 \mu L/min$, the pressure should increase or remain constant at least. If it does not, this indicates possible leakage.

- 7. In this case, find and eliminate the cause for the leak. Possible sources are:
 - Capillary connections
 Inspect them for signs of leakage and perform the test described below this list.
 - *Inline filter in the outlet unit*, especially important with biocompatible pumps Perform the test described below this list.
 - Piston seals
 - ◆ Inspect the piston seals for leakage (→ page 117) and tighten leaking connections, if necessary.
 - If leakage continues, replace the piston seals (\rightarrow page 119).
 - If leakage is observed with new piston seals, operate the pump for at least 2 hours at 350 bar to run in the seals. *Note*: In this case, tightening the pump head screws does not help.
 - Outlet unit seal
 Inspect the seal for signs of leakage and perform the test described below this list.
 - Check valves
 Tighten the check valve screws and perform the test described below this list.
 - ♦ Purge screw

Inspect the purge screw and outlet unit. To make sure that the purge screw is not too tight, perform the test described below this list.

Test

- a) Pressure up the pump. This is the best way to identify a leaking connection.
- b) Allow 5 minutes for the pump to stabilize. This is important as the pressure drops faster during the first 5 minutes, because the seals and other components have to adjust the pressure.
- c) After the stabilization time, monitor the pressure drop.
- d) Tighten the connection a little more tight. The pressure will suddenly increase a little.
- e) Monitor if the pressure drops at the same rate as before. If the pressure drops significantly slower, the connection was leaking.
- 8. Reset the upper pressure limit to the value used before the leakage test.
8 Optimizing the Pump for Special Application

The following pumps have a dynamic mixing chamber (volume: 328 μ L) in the outlet unit: LPG-3400A, LPG-3400AB, DGP-3600A, DGP-3600AB, and HPG-3x00A



Fig. 59: Outlet unit with mixing chamber

| No. | Description |
|-----|--|
| 1 | Outlet unit installation screws |
| 2 | Pressure transducer for the system pressure |
| 3 | Purge valve |
| 4 | Inline filter, consisting of a filter holder and filter frit |
| 5 | Gasket |
| 6 | Stirrer |
| 7 | Mixing chamber (base) |

You may want to optimize the pump for special applications. To do so, the following kits are available:

• Mixing Chamber Extension kits

- to extend the mixing chamber volume by 600 μ L to a total volume of 928 μ L - to extend the mixing chamber volume by 1200 μ L to a total volume of 1528 μ L for the following pumps: LPG-3400A, LPG-3400AB, DGP-3600A, DGP-3600AB, HPG-3200A, and HPG-3400A (\rightarrow page 136)



: As a standard, HPG-3200P pumps are shipped with an extended mixing chamber (total chamber volume = 928μ L).

 Micro Flow kit for pumps LPG-3400A, LPG-3400AB, DGP-3600A, and DGP-3600AB (→ page 138)

8.1 Mixing Chamber Extension

To increase the mixing chamber volume, install the appropriate extension.

| Description | Part No. |
|---|------------------------|
| Mixing chamber extension (to extend the mixing chamber volume by 600 μL) for LPG-3400A, DGP-3600A, HPG-3200A, and HPG-3400A LPG-3400AB and DGP-3600AB The extension kit includes: 1 extension ring (600 μL) 3 screws to attach the extension ring 2 mixing chamber gaskets | 6035.1978 6037.1978 |
| Mixing chamber extension (to extend the mixing chamber volume by 1200 μ L) for LPG-3400A,DGP-3600A, HPG-3200A, HPG-3400A, and HPG-3200P LPG-3400AB and DGP-3600AB The extension kits includes: 1 extension ring (1200 μ L) 3 screws to attach the extension ring 2 mixing chamber gaskets | 6035.1979 6037.1979 |

i Tip: Two kits are required for a DGP-3600A or DGP-3600AB.



Fig. 60: Outlet unit with mixing chamber and optional extension

| No. | Description | No. | Description |
|-----|--|-----|--------------------------|
| 1 | Outlet unit installation screws | 6 | Mixing chamber extension |
| 2 | Pressure transducer (system pressure) | 7 | Gasket |
| 3 | Purge valve | 8 | Stirrer |
| 4 | Inline filter, consisting of a filter holder and filter frit | 9 | Mixing chamber (base) |
| 5 | Gasket | 10 | Outlet unit |

To install the mixing chamber extension

- 1. If necessary, purge the pump to remove toxic solvents.
- 2. Turn off the pump and disconnect the power cord from its source.
- 3. Tilt the front panel upward.
- 4. Disconnect the capillary on the pump outlet.
- 5. Remove the installation screws from the outlet unit (\rightarrow Fig. 60, no. 1), and then carefully remove the outlet unit (\rightarrow Fig. 60, no. 10).



Keep the installation screws in case you wish to remove the extension later.

- 6. Remove the old gasket (\rightarrow Fig. 60, no. 7).
- Place the mixing chamber extension (→ Fig. 60, no. 6) on a flat surface (for example, drip tray). Install the gasket (→ Fig. 60, no. 5) and the outlet unit onto the mixing chamber extension. Insert the long installation screws (shipped with the mixing chamber extension kit) into the openings on the outlet unit.
- Check that the stirrer (→ Fig. 60, no. 8) is in the mixing chamber. Insert the second gasket (→ Fig. 60, no. 7) into the mixing chamber base (→ Fig. 60, no. 9). If the gasket does not adhere to the base, apply some liquid to wet the seal before insertion.
- 9. Hold the outlet unit by the extension ring and insert the assembly into the mixing chamber base. Make sure that the seals (\rightarrow Fig. 60, nos. 5 and 7) remain in their correct positions.
- 10. Install the outlet unit by tightening all installation screws in turn. Tighten the screws just until you suddenly have to use force to turn in the screws. Do not tighten any further to avoid damage to the screws and screwdriver.
- 11. Reconnect the power cord to its source and turn on the pump.
- 12. Purge all channels (\rightarrow page 54).
- 13. Test the pump for leakage (\rightarrow page 133).
- **I** Tip: After you have installed the mixing chamber extension, verify in Chromeleon that the **MixingChamber** property is set to the correct value. If it is not, the diagnostics tests may not provide reliable results.

To uninstall the mixing chamber extension

If you want to remove the extension, follow the steps above. Remove the mixing chamber extension including the ring seal. Reinstall the outlet unit onto the mixing chamber base, using the short installation screws and a new gasket.

8.2 Micro Flow Kit

To reduce the delay volume for a low-pressure pump, install the Micro Flow option.

| Description | Part No. |
|--|------------------------|
| Micro Flow kit for LPG-3400A and DGP-3600A pumps Micro Flow kit for LPG-3400AB and DGP-3600AB pumps The kit contains the following components: 1 L connection 1 Micro flow ring seal 1 capillary from proportioning valve to working head | 6035.1977 6037.1977 |
| Tips: | |
| Two kits are required for a DGP-3600A. LPG-3400M, LPG-3400MB, DGP-3600M, and DGP-3600MB are shipped with the Micro Flow option as standard. | |



Fig. 61: Outlet unit with micro flow option

| No. | Description |
|-----|--|
| 1 | Outlet unit installation screws |
| 2 | Pressure transducer for system pressure |
| 3 | Purge valve |
| 4 | Inline filter, consisting of a filter holder and filter frit |
| 5 | Micro flow ring seal |
| 6 | L connection |
| 7 | Outlet unit |

To install the micro flow option

- 1. If necessary, purge the pump to remove toxic solvents.
- 2. Turn off the pump and disconnect the power cord from its source.
- 3. Tilt the front panel upward.
- 4. Disconnect the capillary from the pump outlet and on the mixing chamber bottom part $(\rightarrow$ Fig. 59, page 135, no. 7).
- 5. Remove the installation screws from the outlet unit (\rightarrow Fig. 61, no. 1), and then carefully remove the outlet unit (\rightarrow Fig. 61, no. 7).
- 6. Remove the mixing chamber base, old gasket, and stirrer (→ Fig. 59, page 135, nos. 5 through 7).
- 7. Install the L connection (\rightarrow Fig. 61, no. 6) to the position of the mixing chamber base so that the capillary connections are in a horizontal position (\rightarrow Fig. 62, page 140).
- 8. Insert the micro flow ring seal (→ Fig. 61, no. 5) into the L connection. If the seal does not adhere to the connection, apply a bit of liquid to wet the seal before insertion.
- 9. Insert the outlet unit into the L connection. Make sure that the seal remains in the correct position.
- 10. Install the outlet unit by tightening all installation screws in turn. Tighten the screws just until you suddenly have to use force to turn in the screws. Do not tighten any further to avoid damage to the screws and screwdriver.
- 11. Reinstall the capillary connections.
- 12. Replace the tubing (ECTFE) from the proportioning valve to the working head
 (→ Fig. 66, page 147 and Fig. 68, page 149) with the PEEK tubing from the Micro Flow option.
 - **I** Tip: Keep the tubing in case you wish to remove the Micro Flow kit later.
- 13. Reconnect the power cord to its source and turn on the pump.
- 14. Purge all channels (\rightarrow page 54).
- 15. Test the pump for leakage (\rightarrow page 133).
- **Tip:** After you have installed the micro flow option, verify in Chromeleon that the **MixingChamber** property is set to the correct value. If it is not, the diagnostics tests may not provide reliable results.

To uninstall the micro flow option

- 1. If necessary, purge the pump to remove toxic solvents.
- 2. Turn off the pump and disconnect the power cord from its source.
- 3. Tilt the front panel upward.
- 4. Disconnect the capillary from the pump outlet and on the L connection.
- 5. Remove the installation screws from the outlet unit (\rightarrow Fig. 61, no. 1), and then carefully remove the outlet unit (\rightarrow Fig. 61, no. 7).
- 6. Remove the L connection (\rightarrow Fig. 61, no. 6) and the old micro flow ring seal.
- 7. Install the mixing chamber base to the position of the L connection so that the capillary connection is in a horizontal position.



Fig. 62: Capillaries connected to the mixing chamber base

- 8. Clean the mixing chamber bottom part and insert the stirrer (\rightarrow Fig. 59, page 135).
- 9. Insert the gasket into the mixing chamber base. If the gasket does not adhere to the base, apply some liquid to wet the seal before insertion.
- 10. Place the outlet unit onto the mixing chamber base. Make sure that the gasket remains in the correct position.
- 11. Install the outlet unit by tightening all installation screws in turn. Tighten the screws just until you suddenly have to use force to turn in the screws. Do not tighten any further to avoid damage to the screws and screwdriver.
- 12. Replace the (PEEK) tubing from the proportioning valve to the working head $(\rightarrow$ Fig. 66, page 147 and Fig. 68, page 149) with the (ECTFE) tubing that was removed during installation of the Micro Flow option.
- 13. Reconnect the capillary to the pump outlet and mixing chamber base.
- 14. Reconnect the power cord to its source and turn on the pump.
- 15. Purge all channels (\rightarrow page 54).
- 16. Test the pump for leakage (\rightarrow page 133).

8.3 Static Mixer

To improve the mixing quality of the two combined solvent streams delivered by a pump HPG-3x00M, install the static mixer.

| Description | Part No. |
|---|-----------|
| Static mixer kit, 50 μL, including: Mixer capillary (0.18 mm I.D.; 180 mm long) + appropriate fittings Static mixer cartridge (50 μL) Static mixer housing (2 pieces) Installation instructions | 6035.0002 |

Install the static mixer as described in the installation instructions shipped with the mixer kit.

Tips: Store the mixer cartridge in a safe place to avoid that the end seals are scratched or damaged.

The mixer cartridge is available also as a separate spare part (part no. 6268.5050).

9 Inside Front Panel Views and Fluid Connections

The sections listed in the table provide a short overview of the inside front panel views and the fluid connections of the UltiMate 3000 system series pumps. In addition, section 9.3.3 describes the special aspects to be considered with a high-pressure gradient system.

| For the | Find information about the | On page |
|------------------------------|---|-------------------|
| Isocratic pump | | |
| ISO-3100A | Inside front panel view Fluid connections | 144 145 |
| Low-pressure gradient pumps | | |
| LPG-3400 | Inside front panel view Fluid connections | 146 147 |
| DGP-3600 | Inside front panel view Fluid connections | 148 149 |
| High-pressure gradient pumps | | |
| HPG-3200 | Inside front panel view Fluid connections | 150 151 |
| HPG-3400 | Inside front panel view Fluid connections | 153 154 |
| Special aspects | Gradient combinations Double Flow mode HPG-3200P semipreparative pump | 156 156 157 |

9.1 Isocratic Pump (IS0-3100A)

9.1.1 Inside Front Panel



Fig. 63: Inside front panel

| No. | Description | Part No. |
|-----|--|--------------------------|
| 1 | Outlet unit with pressure transducer and purge valve $(\rightarrow page 23)$ | |
| 2 | Peristaltic pump | |
| 3 | Leak sensor $(\rightarrow page 25)$ | |
| 4 | Equilibration head | |
| 5 | Working head | |
| 6 | Liquid reservoir of the rear seal wash system | 6030.9501 (2 reservoirs) |

9.1.2 Fluid Connections



Fig. 64: Fluid connections

| No. | Description | Part No. |
|-----|--|------------|
| 1 | Working head | |
| 2 | Capillary from working head to equilibration head | 6030.3015* |
| 3 | Equilibration head | |
| 4 | Capillary from equilibration head to outlet unit | 6035.3011* |
| 5 | Outlet unit with pressure transducer and purge valve $(\rightarrow page 23)$ | |

9.2 Low-Pressure Gradient Pumps

9.2.1 LPG-3400

9.2.1.1 Inside Front Panel



1

2

8 7 Fig. 65: Inside front pan 8 5 4

| No. | Description | Part No. |
|-----|---|-----------------------------|
| 1 | Outlet unit with pressure transducer, purge valve, and inline filter $(\rightarrow page 23)$ It depends on the pump variant whether the outlet unit includes a mixing chamber ($\rightarrow page 24$). | |
| 2 | Peristaltic pump | |
| 3 | Leak sensor (\rightarrow page 25) | |
| 4 | Equilibration head | |
| 5 | Working head | |
| 6 | 4-channel proportioning valve | |
| 7 | Integrated 4-channel vacuum degasser | |
| 8 | Liquid reservoir of the rear seal wash system | 6030.9501 (2 reservoirs) |

9.2.1.2 Fluid Connections

The integrated online degasser is connected to the low-pressure side of the solvent delivery system, that is, the pump inlet.



Fig. 66: Fluid connections (The arrows indicate the direction of flow.)

| No. | Description | Part No. |
|-----|---|--|
| 1 | Integrated online degasser | |
| 2 | Tubing from degasser to proportioning valve | 6035.2540* (set of 4) |
| 3 | 4-channel proportioning valve | |
| 4 | Tubing from proportioning valve to working head LPG -3400A and LPG-3400AB LPG-3400M and LPG-3400MB | 6035.2515* 6035.2514* |
| 5 | Working head | |
| 6 | Capillary from working head to equilibration head LPG-3400A and LPG-3400M LPG-3400AB and LPG-3400MB | 6030.3015* 6037.3015* |
| 7 | Equilibration head | |
| 8 | Capillary from equilibration head to outlet unit LPG-3400A LPG-3400AB LPG-3400M LPG-3400MB | 6035.3011* 6037.3020* 6035.3014* 6037.3011* |
| 9 | Outlet unit with pressure transducer, purge valve, and inline filter $(\rightarrow \text{ page } 23)$. It depends on the pump variant whether the outlet unit includes a mixing chamber $(\rightarrow \text{ page } 24)$. | |

9.2.2 DGP-3600

9.2.2.1 Inside Front Panel



7 6 Fig. 67: Inside front pathel

| No. | Description | Part No. |
|-----|---|-----------------------------|
| 1 | Outlet unit with pressure transducer, purge valve, and inline filter $(\rightarrow page 23)$ It depends on the pump variant whether the outlet unit includes a mixing chamber ($\rightarrow page 24$). | |
| 2 | Working head | |
| 3 | Equilibration head | |
| 4 | Peristaltic pump | |
| 5 | Leak sensor $(\rightarrow page 25)$ | |
| 6 | 3-channel proportioning valve | |
| 7 | Liquid reservoir of the rear seal wash system | 6030.9501 (2 reservoirs) |



9.2.2.2 Fluid Connections

Fig. 68: Fluid connections (The arrows indicate the direction of flow.)

| No. | Description | Part No. |
|-----|---|--|
| | SRD-3600 Solvent Rack with analytical 6-channel vacuum degasser | 5035.9230 |
| 1 | Solvent lines from degasser to proportioning valve | 6030.2547* (set of 3) |
| 2 | 3-channel proportioning valve | |
| 3 | Tubing from proportioning valve to working head DGP-3600A and DGP-3600AB DGP-3600M and DGP-3600MB | 6035.2515* 6035.2514* |
| 4 | Working head | |
| 5 | Capillary from working head to equilibration head DGP-3600A and DGP-3600M DGP-3600AB and DGP-3600MB | 6030.3015* 6037.3015* |
| 6 | Equilibration head | |
| 7 | Capillary from equilibration head to outlet unit DGP-3600A DGP-3600AB DGP-3600M DGP-3600MB | 6035.3011* 6037.3020* 6035.3014* 6037.3011* |
| 8 | Outlet unit with pressure transducer, purge valve, and inline filter $(\rightarrow \text{ page } 23)$ It depends on the pump variant whether the outlet unit includes a mixing chamber (\rightarrow page 24). | |

9.3 High-Pressure Gradient Pumps

9.3.1 HPG-3200

9.3.1.1 Inside Front Panel



Fig. 69: Inside front panel (here: HPG-3200 with mixing chamber in the outlet unit)

| No. | Description | Part No. |
|-----|---|-----------------------------|
| 1 | Tee insert (HPG-3200A and HPG-3200P only) | 6035.1980 |
| 2 | Working head with pressure transducer | |
| 3 | Equilibration head | |
| 4 | Outlet unit with pressure transducer, purge valve, mixing chamber, and inline filter (\rightarrow page 23) | |
| 5 | Peristaltic pump | |
| 6 | Leak sensor (\rightarrow page 25) | |
| 7 | Liquid reservoir of the rear seal wash system | 6030.9501 (2 reservoirs) |

An HPG-3200M has no mixing chamber. Thus, the equilibration head of the left pump block is connected directly to the outlet unit (\rightarrow Fig. 71, page 152). Due to limited space, it may be difficult to install the capillary from the left equilibration head to the outlet unit. To facilitate the installation, follow the steps on page 129.

A static mixer is available as an option for the HPG-3200M pump (\rightarrow page 141). The mixer improves the mixing quality of the 2 combined solvent streams delivered by the pump, and thus guarantees a smoother baseline.

9.3.1.2 Fluid Connections



Fig. 70: Fluid connection (HPG-3200A with mixing chamber in the outlet unit) The arrows indicate the direction of flow. With an HPG-3200P, the capillaries run slightly different.

| No. | Description | Part No. |
|-----|---|--|
| | SRD-3200 Solvent Rack with analytical 2-channel vacuum degasser (for use with HPG-3200A only) | 5035.9250 |
| 1 | Solvent lines from analytical degasser to working head (for use with HPG-3200A only) | 6035.2530* (set of 2) |
| 2 | Working head with pressure transducer | |
| 3 | Capillary from working head to equilibration head HPG-3200A HPG-3200P | 6030.3015* 5030.3025* |
| 4 | Equilibration head | |
| 5 | Capillary from left equilibration head to Tee insert HPG-3200A HPG-3200P | Included in: 6035.1967* 6035.1969* |
| 6 | Capillary from right equilibration head to Tee insert HPG-3200A HPG-3200P | Included in: 6035.1967* 6035.1969* |
| 7 | Tee insert | 6035.1980 |
| 8 | Capillary from Tee insert to outlet unit HPG-3200A HPG-3200P | Included in: 6035.1967* 6035.1969* |
| 9 | Outlet unit with pressure transducer, purge valve, mixing chamber, and inline filter (\rightarrow page 23) | |



Fig. 71: Fluid connections (HPG-3200M) The arrows indicate the direction of flow.

| No. | Description | Part No. |
|-----|--|----------------------------|
| | SRD-3200 Solvent Rack with analytical 2-channel vacuum degasser | 5035.9250 |
| 1 | Solvent lines from degasser to working head | 6035.2530* (set of 2) |
| 2 | Working head with pressure transducer | |
| 3 | Capillary from working head to equilibration head | 6030.3015* |
| 4 | Equilibration head | |
| 5 | Capillary from left equilibration head to outlet unit | Included in: 6035.1968* |
| 6 | Capillary from right equilibration head to outlet unit | Included in: 6035.1968* |
| 7 | Outlet unit with pressure transducer, purge valve, and inline filter $(\rightarrow page 23)$ | |

9.3.2 HPG-3400



Fig. 72: Inside front panel (here: HPG-3400A with mixing chamber in the outlet unit)

| No. | Description | Part No. |
|-----|---|-----------------------------|
| 1 | Tee insert (HPG-3400A only) | 6035.1980 |
| 2 | Working head with pressure transducer | |
| 3 | Equilibration head | |
| 4 | Outlet unit with pressure transducer, purge valve, mixing chamber, and inline filter (\rightarrow page 23) | |
| 5 | Peristaltic pump | |
| 6 | Leak sensor (\rightarrow page 25) | |
| 7 | Solvent Selector "2 from 4" | |
| 8 | Liquid reservoir of the rear seal wash system | 6030.9501 (2 reservoirs) |

An HPG-3400M has no mixing chamber. Thus, the equilibration head of the left pump block is connected directly to the outlet unit (\rightarrow Fig. 74, page 155). Due to limited space, it may be difficult to install this capillary from the left equilibration head to the outlet unit. To facilitate the installation, follow the steps on page 129.

A static mixer is available as an option for the HPG-3200M pump (\rightarrow page 141). The mixer improves the mixing quality of the 2 combined solvent streams delivered by the pump, and thus guarantees a smoother baseline.

9.3.2.2 Fluid Connections



Fig. 73: Fluid connections (HPG-3400A with mixing chamber in the outlet unit) The arrows indicate the direction of flow.

| No. | Description | Part No. |
|-----|---|----------------------------|
| | SRD-3400 Solvent Rack with analytical 4-channel vacuum degasser | 5035.9245 |
| 1 | Solvent lines from degasser to solvent selector | 6035. 2532* (set of 4) |
| 2 | Solvent selector "2 from 4" | |
| 3 | Tubing from solvent selector to working head | 6035.2515* |
| 4 | Working head with pressure transducer | |
| 5 | Capillary from working head to equilibration head | 6030.3015* |
| 6 | Equilibration head | |
| 7 | Capillary from left equilibration head to Tee insert | Included in: 6035.1967* |
| 8 | Capillary from right equilibration head to Tee insert | Included in: 6035.1967* |
| 9 | Tee insert | 6035.1980 |
| 10 | Capillary from Tee insert to outlet unit | Included in: 6035.1967* |
| 11 | Outlet unit with pressure transducer, purge valve, mixing chamber, and inline filter (\rightarrow page 23) | |



Fig. 74: Fluid connections (HPG-3400M) The arrows indicate the direction of flow.

| No. | Description | Part No. |
|-----|--|--------------------------|
| | SRD-3400 Solvent Rack with analytical 4-channel vacuum degasser | 5035.9245 |
| 1 | Solvent lines from degasser to solvent selector | 6035.2532* (set of 4) |
| 2 | Solvent selector "2 from 4" | |
| 3 | Tubing from solvent selector to working head | 6035.2515* |
| 4 | Working head with pressure transducer | |
| 5 | Capillary from working head to equilibration head | 6030.3015* |
| 6 | Equilibration head | |
| 7 | Capillary from left equilibration head to outlet unit | Included in 6035.1968* |
| 8 | Capillary from right equilibration head to outlet unit | Included in 6035.1968* |
| 9 | Outlet unit with pressure transducer, purge valve, and inline filter $(\rightarrow page 23)$ | |

9.3.3 Special Aspects of a High-Pressure Gradient System

The **HPG-3x00** high-pressure gradient pumps operate with two pump blocks that are integrated in a single enclosure. Operation of the high-pressure gradient pump is identical to the low-pressure gradient pump (LPG). However, consider the information in the following sections.

9.3.3.1 Supported Gradient Combinations

For an HPG-3400 pump with "2 from 4" solvent selector, channels A and C are connected to pump block 1 (left), while channels B and D are connected to pump block 2 (right). The solvent selector supports the following combinations of binary high-pressure gradients:

| Pump block 1 (left) | Pump block 2 (right) |
|---------------------|----------------------|
| Solvent A | Solvent B |
| Solvent A | Solvent D |
| Solvent C | Solvent B |
| Solvent C | Solvent D |

In a specific time step, each pump block delivers only one of the two channels (100%). It is not possible to specify a mixing ratio for the two channels connected to the same pump block (A and C; B and D).

If you enter a portion for solvent C while solvent A is already set, the portion for solvent A is automatically reduced to 0%. The same applies to channels B and D on the second pump block. For example, if you enter a portion for solvent D (D = 20%), while the portions for solvents B and C are already set (B = 40%, C = 60%), solvent C is automatically increased to 80%. The same applies to all other channels.

9.3.3.2 Double Flow Mode

HPG models support the double flow mode. This mode enables maximum flow rates of 20 mL/min (analytical and micro pumps) or 100 mL/min (semipreparative pump). The double flow mode is available for each solvent combination and composition supported by the HPG-pump. The pump automatically calculates the flow rate maximum considering that none of the two pump blocks delivers more than 10 mL/min (analytical version) or 50 mL/min (semipreparative pump).

Example A:

Both pump blocks of an analytical pump deliver 50%. Each block then delivers 10 mL/min thus, reaching the flow rate maximum of 20 mL/min.

Example B:

Pump block A of an analytical pump delivers 60% and block B delivers 40%. Block A then delivers 10 mL/min while block B delivers correspondingly less, that is, 6.667 mL/min. In this case, the total maximum flow rate is 16.667 mL/min.

9.3.4 HPG-3200P Semipreparative Pump

The HPG-3200P is the semipreparative pump in the UltiMate 3000 system series. Please observe the following when operating this pump:

- Only use the solvent lines shipped with the pump (part no. 6250.4301). These lines have an inner diameter of 3.0 mm.
- Only use the filter frits from the accessories kit (also included in part no. 6035.1964).
- Do not prolong the solvent lines.
- Especially for applications with high flow rates, make sure that the solvent reservoirs are located near the pump; they should be on the same level or higher. This is to avoid formation of air bubbles in the reservoirs.
- Thermo Fisher Scientific recommends using a preparative degasser. Do not connect an analytic degasser to the semipreparative pump. Else, degas the solvent, for example, in an ultrasonic bath.

10 Technical Information

Technical information: November 201. All technical specifications are subject to change without notice.

| | ISO-3100A | LPG-3400A/ LPG-3400AB | DGP-3600A DGP-3600AB | HPG-3200A/ HPG-3400A | HPG-3200P | |
|--|---|---|---|----------------------------------|---|--|
| Flow rate range Recommended Settable | μL/min 50–10,000 1-10,000 | μL/min 200–10,000 1-10,000 | μL/min 200–10,000 1-10,000 | μL/min 100–10,000 1-10,000 | mL/min 0.5–50 0.001-50 | |
| Flow rate accuracy | | at l | ±0.1% mL/min | | ±0,1% | |
| Flow rate precision | < | 0.05% RSD or <0.01 | min SD, whichever is | greater | <0.1% RSD at 3 mL/min | |
| Pressure range | | 2-50 M | Pa (7250 psi) | | 1-10 MPa (1450 psi) (short term: 15 MPa-2100 psi) | |
| Pressure ripple | | Typical <0.2 MPa or | <1%, whichever is gro | eater | Typical <0.2 MPa or <1% whichever is greater | |
| Proportioning accuracy | | ±0.5% of fullscale | ±0.5% of fullscale | ±0.2% of fullscale | ±0.2% of fullscale | |
| Proportioning precision | | <0.2% SD at 2 mL/min | <0.2% SD at 2 mL/min | <0.2% SD at 2 mL/min | <0.2% SD at 5 mL/min | |
| No of eluent lines | nt lines 1 | | 6 (2 x 3) | 2 | 2 | |
| Grad. delay volume | | 690 μL or LPG-3400A: 360 μL with Micro Flow kit | 690 μL or DGP-3600A: 360 μL with Micro Flow kit | 375 μL | 1035 μL | |
| Solvent degassing | External (optional) | Built-in 4-channel degasser (channel volume: 670 µL) | External (optional) | External (optional) | | |
| Weight | 16.2 kg (35.2 lb) 17.9 kg (39.4 lb) 21.5 kg (47.3 lb) 20.8 kg (45.8 lb) 20.8 kg | | | | 20.8 kg (45.8 lb) | |
| Dimensions | (h x w x d): 19 x 4 | ¹ 2 x 51 cm (7.5 x 16.5 | x 20 in.) | | | |
| PC connection | All functions controllable via USB 1.1; integrated USB hub with three free USB ports | | | | | |
| I/O interfaces | 3 digital inputs, 4 15-pin D-Sub con | 4 programmable relays, 1 analog output for pressure monitoring inector for Solvent Rack/degasser connection | | | | |
| Safety features | Leak sensor Active rear seal w reservoir | eak sensor ctive rear seal washing with monitoring of the piston seal tightness and the liquid level of the liquid servoir | | | | |
| User input/ display | LCD indicating sy Standby button; 3 4 function keys fo | system parameters 3 LEDs (Power, Connected, and Status) for status monitoring for operation during initial installation and maintenance | | | | |
| GLP | In Chromeleon: Full su monitoring. All system | | Il support of automatic equipment qualification (AutoQ [™]) and System Wellness stem parameters are logged in the Chromeleon Audit Trail. | | | |
| Wetted parts | Stainless steel or titanium (for biocompatible pumps), sapphire, ruby, UHMW polyethylene, PCTFE, PTFE, PEEK™, zirconium oxide, aluminum oxide, FEP, ECTFE | | | | | |
| Power requirements | 150 VA (100-120V, 60 Hz; 200-240V, 50 Hz) | | | | | |
| Emission sound pressure level: | < 70 dB(A) in 1-m-distance | | | | | |
| Environmental conditions | Range of use: Indoor use; Temperature: 10 °C to 35 °C; Air humidity:80% relative humidity, non-condensing; Overvoltage category: II; Pollution degree: 2 | | | | | |

| | HPG-3200M/ HPG-3400M | LPG-3400M LPG-3400MB | DGP-3600M/ DGP-3600MB | LPG-3400M(B) + FLM-3x00(B) | DGP-3600M(B) + FLM-3x00(B) | |
|--|--|---|---|---|---|--|
| Flow rate range Recommended (Settable) | µL/min 20–10,000 (1-10,000) | μL/min 10–2,500 (1-2,500) | μL/min 10–2,500 (1-2,500) | Nano: 50–1,000 nL/min Cap: 0.5-10 μL/min Micro: 10-160 μL/min | <i>Nano:</i> 50–1,000 nL/min <i>Cap:</i> 0.5-10 μL/min <i>Micro:</i> 10-160 μL/min | |
| Flow rate accuracy | ±0.1% at 1 mL/min | ±0.5% at 200 μL/min (typical ±0.25% at 200 μL/min) | ±0.5% at 200 μL/min (typical ±0.25% at 200 μL/min) | Nano: ±3% at 300 nl/min Cap: ±1.5% at 4 μL/min Mikro: ±1% at 50 μL/min | Nano: ±3% at 300 nl/min Cap: ±1.5% at 4 μL/min Mikro: ±1% at 50 μL/min | |
| Flow rate precision | <0.05% RSD or | <0.01 min SD, whichev | ver is greater | <0.1% RSD | <0.1% RSD | |
| Pressure range | | 2-50 MPa (7250 psi) | | 1-50 MP | a (7250 psi) | |
| Pressure ripple | Tyj | pical <0.2 MPa or <1% whichever is greater | | | | |
| Proportioning accuracy | | $\pm 0.2\%$ of fullscale | | $\pm 2\%$ of fulls | cale on column | |
| Proportioning precision | <0.2% SD at 2 mL/min | <0.5% SD at 200 µL/min | <0.5% SD at 200 µL/min | <0.5% SD on column | <0.5% SD on column | |
| No. of eluent lines | 2 | 4 | 6 (2 x 3) | 4 | 6 (2 x 3) | |
| Grad. delay volume | 35 µL | 230 µL | 230 µL | Effektiv: Nano: 0.52 μL Cap: 3.3 μL Micro: 50 μL | Effektiv: <i>Nano:</i> 0.52 μL <i>Cap:</i> 3.3 μL <i>Micro:</i> 50 μL | |
| Solvent degassing | External (optional) | Built-in 4-channel degasser (channel volume: 670 µL) | External (optional) | Built-in 4-channel degasser (channel volume: 670 µL) | External (optional) | |
| Weight | 20.8 kg (45.8 lb) | 17.9 kg (39.4 lb) | 21.5 kg (47.3 lb) | <i>LPG:</i> 17.9 kg (39.4 <i>lb</i>) <i>FLM:</i> 17 kg (37.4 lb) | DGP: 21.5 kg (47.3 lb) FLM: 17 kg (37.4 lb) | |
| Dimensions | (h x w x d): 19 x 42 x 5 | 51 cm (7.5 x 16.5 x 20 ir | 1.) | | | |
| PC connection | All functions controllal | ble via USB 1.1; integra | ted USB hub with th | nree free USB ports | | |
| I/O interfaces | 3 digital inputs, 4 programmable relays, 1 analog output for pressure monitoring 15-pin D-Sub connector for Solvent Rack/degasser connection | | | | | |
| Safety features | Leak sensor Active rear seal washing with monitoring of the piston seal tightness and the liquid level of the liquid reservoir | | | | | |
| User input/ display | LCD indicating system parameters, Standby button, 3 LEDs (Power, Connected, and Status) for statu monitoring, 4 function keys for operation during initial installation and maintenance | | tus) for status | | | |
| GLP | In Chromeleon: Full support of automatic equipment qualification (AutoQ TM) and System Wellness monitoring. All system parameters are logged in the Chromeleon Audit Trail. | | | | | |
| Wetted parts | Stainless steel or titanium (for biocompatible pumps), sapphire, ruby, UHMW polyethylene, PCTFE, PTFE, PEEK™, zirconium oxide, aluminum oxide, FEP, ECTFE | | | | | |

| | HPG-3200M/ HPG-3400M | LPG-3400M LPG-3400MB | DGP-3600M/ DGP-3600MB | LPG-3400M(B) + FLM-3x00(B) | DGP-3600M(B) + FLM-3x00(B) |
|--|---|-------------------------|--------------------------|-------------------------------|-------------------------------|
| Power requirements | 150 VA (100-120V, 60 Hz; 200-240V, 50 Hz) | | | | |
| Emission sound pressure level: | < 70 dB(A) in 1-m-distance | | | | |
| Environmental conditionsRange of use: Indoor use; Temperature: 10 °C to 35 °C; Air humidity:80% relative humidity, non-condensing; Overvoltage category: II; Pollution degree: 2 | | | dity, | | |

11 Accessories, Spare Parts, and Consumables

Accessories, spare parts, and consumables for the pump are always maintained at the latest technical standard. Therefore, part numbers are subject to alteration. However, updated parts will always be compatible with the parts they replace.

11.1 Standard Accessories

The following accessories are shipped with the pump. (Note: The list is subject to change without notice.) Some parts listed in the following tables are included in one of the spare part kits. For information about these kits, see section 11.3 (\rightarrow page 169).

The part number always refers to the packing unit. Unless otherwise stated, the packing unit is 1 unit. For more information, contact the Thermo Fisher Scientific sales organization for Dionex HPLC Products

| Description | Part No. | Quantity in the accessories kit |
|---|--|-----------------------------------|
| Accessory kit for ISO-3100A, LPG-3400A, LPG-3400M, DGP-3600A, DGP-3600M, HPG-3x00A, and HPG-3x00M pumps, including | | |
| Filter frit for inline filter (stainless steel, porosity: $0.5 \ \mu m$) | Included in 6000.0045 | 2 filter frits |
| Filter frit for inline filter (stainless steel, porosity: 10 µm) | Included in 6268.0031 | 2 filter frits |
| Solvent filter, including: Filter holder (top and bottom parts) and filter frit (stainless steel, 10 µm) | Included in 6268.0115 Included in 6268.0110 | 1 |
| Tool kit for analytical UltiMate 3000 pumps, with 1 off: Double open-end wrench $(1/4 \times 5/16)$, double open-end wrench (11×13) , Allen wrench (size 3) | Included in 6007.9302 | 1 |
| Capillary (1/16", 0.25 mm ID) with appropriate fittings | 6020.9110 | 1 |
| Menu pen | 6300.0100 | 1 |
| Fuse, 2A, TT (5 x 20 mm) | Included in 6030.9003 | 2 |
| Silicone tubing (OD x ID: 2.80 x 1.30 mm) | Included in 6000.0010 | 3 m |
| Components for UltiMate 3000 system drainage: Cable clips (self-adhesive) Y piece T piece Elbow Connecting tube Drain tubing Installation instructions | Included in 6040.0005 | 4 5 4 12 1 6m 1 |

| Description | Part No. | Quantity in the accessories kit |
|---|-----------------------|---------------------------------|
| Set of PEEK fittings (for tubing O.D. 3.0 x I.D. 1.5 mm), including 5 fittings (1/8" PEEK) 5 locking rings (PEEK) 5 support flanges (PEEK) 2m solvent supply line (analytical) | 6025.2511 | 1 |
| Piston support hemisphere | 6269.0040 | 2 |
| Solvent reservoir (250 mL) with bottle cap | 2270.0026 | 1 |
| Seal wash reservoir for unidirectional seal washing | | 1 |
| Plastic syringe (12 mL) | Included in 6000.0010 | 1 syringe |
| Extractor for rear seal wash body | 6000.0046 | 1 |
| HPLC Troubleshooting Guide Poster | 6040.0050 | 1 |
| USB cable, type A to type B, USB 2.0, 1 m | 6035.9035 | 1 |
| USB cable, type A to type B, USB 2.0, 5 m | 6911.0002 | 1 |

| Description | Part No. | Quantity in the accessories kit |
|---|-----------------------|-----------------------------------|
| Accessory kit for LPG-3400AB, LPG-3400MB, DGP-3600AB and DGP-3600MB pumps, including: | | |
| Filter frit for inline filter (titanium, porosity: 2 µm) | Included in 6268.0036 | 2 filter frits |
| Filter frit for inline filter (titanium, porosity: 10 µm) | Included in 6268.0032 | 2 filter frits |
| Tool kit for analytical UltiMate 3000 pumps, with 1 off: Double open-end wrench $(1/4 \times 5/16)$, double open-end wrench (11×13) , Allen wrench (size 3) | Included in 6007.9302 | 1 |
| Capillary (PEEK, OD x ID: 1/16" x 0.25 mm) | 6251.6001 | 2 m |
| Fitting screw (PEEK, 1/16", 15 mm) | 6266.0024 | 1 |
| Menu pen | 6300.0100 | 1 |
| Fuse, 2A, TT (5 x 20 mm) | Included in 6030.9003 | 2 |
| Silicone tubing (OD x ID: 2.80 x 1.30 mm) | Included in 6000.0010 | 3 m |
| Components for UltiMate 3000 system drainage: Cable clips (self-adhesive) Y piece T piece Elbow Connecting tube Drain tubing Installation instructions | Included in 6040.0005 | 4 5 4 12 1 6m 1 |

| Description | Part No. | Quantity in the accessories kit |
|---|--|---------------------------------|
| Set of PEEK fittings (for tubing O.D. 3.0 x I.D. 1.5 mm), including 5 fittings (1/8" PEEK) 5 locking rings (PEEK) 5 support flanges (PEEK) 2m solvent supply line (analytical) | 6025.2511 | 1 |
| Solvent line filter, including: Filter holder (top and bottom parts) Filter frit (Titanium, porosity: 10 μm) | Included in 6268.0115 Included in 6268.0111 | 1 6 |
| Piston support hemisphere | 6269.0040 | 2 |
| Solvent reservoir (250 mL) with bottle cap | 2270.0026 | 1 |
| Seal wash reservoir for unidirectional seal washing | | 1 |
| Plastic syringe (12 mL) | Included in 6000.0010 | 1 syringe |
| Extractor for rear seal wash body | 6000.0046 | 1 |
| HPLC Troubleshooting Guide Poster | 6040.0050 | 1 |
| USB cable, type A to type B, USB 2.0, 1 m | 6035.9035 | 1 |
| USB cable, type A to type B, USB 2.0, 5 m | 6911.0002 | 1 |

| Description | Part No. | Quantity in the accessories kit |
|---|-----------------------|-----------------------------------|
| Accessory kit for HPG-3200P pumps, including: | | |
| Filter frit for inline filter (stainless steel, porosity: $10 \ \mu m$) | Included in 6268.0031 | 2 filter frits |
| Capillary (1/16", 0.70 mm ID) with appropriate fittings | 6005.9100 | 1 |
| Tool kit for semipreparative pump, with 1 off: Double open-end wrench $(1/4 \times 5/16)$, double open-end wrench (11×13) , Allen wrench (size 3), slotted screwdriver (for 8mm pistons), installation tool for cap seal | 6007.9303 | 1 |
| Menu pen | 6300.0100 | 1 |
| Fuse, 2A, TT (5 x 20 mm) | Included in 6030.9003 | 2 |
| Silicone tubing (OD x ID: 2.80 x 1.30 mm) | Included in 6000.0010 | 3 m |
| Components for UltiMate 3000 system drainage: Cable clips (self-adhesive) Y piece T piece Elbow Connecting tube Drain tubing Installation instructions | Included in 6040.0005 | 4 5 4 12 1 6m 1 |

| Description | Part No. | Quantity in the accessories kit |
|--|-----------------------|---------------------------------|
| Solvent line filter (preparative) with Tubing adaptor (ID 3.0 to ¹ / ₄ -28") | 6035.1964 | 2 |
| Solvent line (3.0 mm ID) | 6250.4301 | 2 |
| Solvent reservoir (250 mL) with bottle cap | 2270.0026 | 1 |
| Seal wash reservoir for unidirectional seal washing | | 1 |
| Plastic syringe (12 mL) | Included in 6000.0010 | 1 syringe |
| Extractor for rear seal wash body | 6000.0046 | 1 |
| HPLC Troubleshooting Guide Poster | 6040.0050 | 1 |
| USB cable, type A to type B, USB 2.0, 1 m | 6035.9035 | 1 |
| USB cable, type A to type B, USB 2.0, 5 m | 6911.0002 | 1 |

11.2 Optional Accessories

| Accessory | Description | Part No. |
|---|---|------------------------|
| Adaptor and Blind Nut kit | <i>For biocompatible pumps</i> The kit is required in addition to the Diagnostics Tool kit if you want to run the diagnostics tests for the biocompatible pumps of the UltiMate 3000 system. The kit includes 1 adaptor capillary (0.3 m; 1.58 mm x 0.5 mm AD x ID) incl. appropriate fitting connections, 2 blind nuts (FS-8), 1 coupling piece (1(16"), and 2 ferrules (1/16" PEEK) | 6037.3012 |
| Diagnostics Tool kit | <i>For all pumps except HPG-3200P</i> The kit is required, for example, when performing certain diagnostics tests with Chromeleon. The kit includes a backpressure capillary (15 m, 0.18 mm ID) including appropriate fittings and two blind nuts (FS-8). | 6035.3000 |
| Manual injection valve | Including installation kit Analytical/micro Semipreparative | 5035.0600 5035.0601 |
| Manual injection valve (installation kit) | The kit does <i>not</i> include the manual injection valve. | 6035.0610 |
| Micro Flow kit | Reduces the delay volume For LPG-3400A and DGP-3600A pumps For LPG-3400AB and DGP-3600AB pumps | 6035.1977 6037.1977 |
| Mixing Chamber Extension kit | <i>For LPG-3400A, DGP-3600A, HPG-3200A, and HPG-3400A pumps</i> Extends the mixing chamber volume by 600 μL Extends the mixing chamber volume by 1200 μL | 6035.1978 6035.1979 |
| | <i>For HPG-3200P pumps</i> Extends the mixing chamber volume by 1200 μL | 6035.1979 |
| | For LPG-3400AB and DGP-3600AB pumps Extends the mixing chamber volume by 600 μL Extends the mixing chamber volume by 1200 μL | 6037.1978 6037.1979 |
| Syringe for use with manual injection valve | 100 μ L syringe for use with analytical/micro valve 5 mL syringe for use with semipreparative valve | 6035.0665 6035.0670 |
| SR-3000 | Solvent Rack without vacuum degasser For use with an LPG-3400A(B) or LPG-3400M(B) pump. | 5035.9200 |
| SRD-3200 | Solvent Rack with analytical 2-channel vacuum degasser For use with an ISO-3100Aor HPG-3200A or M pump. | 5035.9250 |
| SRD-3400 | Solvent Rack with analytical 3-channel vacuum degasser For use with an HPG-3400A or M pump or with two HPG-3200A or M pumps in a two-stack system | 5035.9245 |

| Accessory | Description | Part No. |
|------------------|---|------------------------|
| SRD-3600 | Solvent Rack with analytical 6-channel vacuum degasser For use with a DGP-3600 (A(B) or M(B)) pump or with one HPG-3400A or M and one HPG-3200A or M pump in a two- stack system | 5035.9230 |
| Static Mixer kit | The static mixer improves the mixing quality of the 2 combined solvent streams delivered by an HPG-3x00M pump, and thus guarantees a smoother baseline that may be required by special applications. The kit includes the mixer housing (2 part), mixer cartridge, and capillary including the appropriate fittings. | 6035.0002 6268.5050 |
| | The mixer cartridge is available also as a separate part. | 0200.0000 |

11.3 Consumables and Spare Parts

The part number always refers to the packing unit. Unless otherwise stated, the packing unit is 1 unit. For more information, contact the Thermo Fisher Scientific sales organization for Dionex HPLC Products

| Description | Part No. |
|---|-----------|
| Adaptor and Blind Nut kit (for biocompatible pumps; \rightarrow section 11.2, page 167) | 6037.3012 |
| Blind nut (FS-8; for all pumps except biocompatible pumps) | 6000.0044 |
| Blind nut (FS-8 biocompatible; for biocompatible pumps) | 6000.0144 |
| Capillary (1/16", 0.25 mm ID) with appropriate fittings (for ISO-3100A, LPG-3400A, LPG-3400M, DGP-3600A, DGP-3600M, HPG-3x00A, and HPG-3x00M pumps) | 6020.9110 |
| Capillary (1/16", 0.70 mm ID) with appropriate fittings (for HPG-3200P) | 6005.9100 |
| Capillary (PEEK, OD x ID: 1/16" x 0.25 mm) (for LPG-3400AB, LPG-3400MB, DGP-3600AB, and DGP-3600MB) | 6251.6001 |
| Capillary (long) from loading pump to autosampler switching valve, standard devices (130 µm ID x 75 cm, PEEK, including appropriate fittings | 6720.0032 |
| Capillary (long) from loading pump to autosampler switching valve, biocompatible devices (130 μ m ID x 75 cm, PEEK, including appropriate fittings for biocompatible pumps) | 6721.0032 |
| Capillary (short) from loading pump and to autosampler switching valve (standard devices), (130 µm ID x 60 cm, PEEK, including appropriate fittings) | 6720.0031 |
| Capillary (short) from loading pump and to autosampler switching valve (biocompatible devices), (130 μ m ID x 60 cm, PEEK, including appropriate fittings | 6721.0031 |
| Capillary (long) from pump to flow manager (standard devices) for 2D LC Comprehensive applications, including the appropriate fittings and ferrules | 6035.2556 |
| Capillary (long) from pump to flow manager (biocompatible devices) for 2D LC Comprehensive applications, including the appropriate fittings and ferrules | 6037.2556 |
| Capillary (short) from pump to flow manager, standard devices, for 2D LC Comprehensive applications, including the appropriate fittings and ferrules | 6035.2554 |
| Capillary (short) from pump to flow manager, biocompatible devices, for 2D LC Comprehensive applications, including the appropriate fittings and ferrules | 6037.2554 |
| Capillary (long) from pump to flow manager (standard devices), including the appropriate fittings and ferrules | 6035.2550 |
| Capillary (long) from pump to flow manager (biocompatible devices), including the appropriate fittings and ferrules | 6037.2550 |
| Capillary (short) from pump to flow manager (standard devices), including the appropriate fittings and ferrules | 6035.2553 |
| Capillary (short) from pump to flow manager (biocompatible devices), including the appropriate fittings and ferrules | 6037.2553 |

| Description | Part No. |
|---|-----------|
| Capillary from working head to equilibration head (IS0-3100A, LPG-3400A(B), LPG-3400M, DGP-3600A(B), DGP-3600M, HPG-3x00A, and HPG-3x00M), including the appropriate fittings and ferrules | 6030.3015 |
| Capillary from working head to equilibration head (LPG-3400AB and MB and DGP-3600AB and MB), including the appropriate fittings and ferrules | 6037.3015 |
| Capillary from working head to equilibration head (HPG-3200P), including the appropriate fittings and ferrules | 5030.3025 |
| Capillary from equilibration head to outlet unit (ISO-3100A, LPG-3400A, and DGP-3600A) | 6035.3011 |
| Capillary from equilibration head to outlet unit (LPG-3400M and DGP-3600M) | 6035.3014 |
| Capillary from equilibration head to outlet unit (LPG-3400MB and DGP-3600MB) | 6037.3011 |
| Capillary from equilibration head to outlet unit (LPG-3400AB and DGP-3600AB) | 6037.3020 |
| Capillary kit (HPG-3x00A), including: Capillary from left equilibration head to Tee insert, Capillary from right equilibration head to Tee insert, Capillary from Tee insert to outlet of the nit outlet including the appropriate fittings and ferrules | 6035.1967 |
| Capillary kit (HPG-3x00M), including: Capillary from left equilibration head to outlet unit, Capillary from right equilibration head to outlet unit including the appropriate fittings and ferrules | 6035.1968 |
| Capillary kit (HPG-3200P), including: Capillary from left equilibration head to Tee insert, Capillary from right equilibration head to Tee insert, Capillary from Tee insert to outlet of the outlet unit including the appropriate fittings and ferrules | 6035.1969 |
| Check valve cartridge (sapphire), (to be used with all pump types, including biocompatible pump versions) | 6035.2300 |
| Check valve nut kit for double check valve for ISO-3100A, LPG-3400A or M, DGP-3600A or M, and HPG-3x00A or M, including: inlet check valve nut and outlet check valve nut | 6035.1965 |
| Check valve nut kit for double check valve for LPG-3400AB and MB and DGP-360AB and MB, including inlet check valve nut and outlet check valve nut | 6037.1964 |
| Check valve nut kit for double check valve for HPG-3200P, including: inlet check valve nut and outlet check valve nut | 6035.1966 |
| Diagnostics Tool kit (for all pumps except HPG-3200P; \rightarrow section 11.2, page 167) | 6035.3000 |
| Drain kit for UltiMate 3000 systems The kit includes all required components and detailed installation instructions. | 6040.0005 |
| Filter frit for inline filter (outlet unit) The kit includes 10 titanium frits (porosity:2 μm). | 6268.0036 |
| Filter frit for inline filter (outlet unit) The kit includes 10 stainless steel frits (porosity: 0.5 μm). | 6000.0045 |
| Description | Part No. |
|---|-----------|
| Filter frit for inline filter (outlet unit) The kit includes 10 stainless steel frits (porosity: 10 µm). | 6268.0031 |
| Filter frit for inline filter (outlet unit) The kit includes 10 titanium frits (porosity: 10 µm). | 6268.0032 |
| Filter frit for solvent line filter The kit includes 10 titanium filter frits (porosity: 10µm). | 6268.0111 |
| Filter frit for solvent line filter The kit includes 10stainless steel filter frits (porosity: 10 μm). | 6268.0110 |
| Filter frit kit (preparative) for HPG-3200P The kit includes 2 filter frits (stainless steel) and 2 tubing adaptors (ID 3.0 to ¹ / ₄ -28") | 6035.1964 |
| Filter holder (PEEK) for solvent line filter, including: 6 filter holder bottom parts and 6 filter holder top parts | 6268.0115 |
| Filter holder (outlet unit), stainless steel pump | 6030.2103 |
| Filter holder (outlet unit), biocompatible pump | 6030.2112 |
| Filter removal tool (1/4"), for the filter in the filter holder on the outlet unit | 6035.3051 |
| Fitting screw (PEEK, 1/16", 15 mm) (for LPG-3400AB, LPG-3400MB, DGP-3600AB, and DGP-3600MB) | 6266.0024 |
| Fitting set for tubing (OD 3.0 x ID 1.5 mm), including: 2 solvent lines (OD 3.0 x ID 1.5 mm), 5 fittings (1/8", PEEK), 5 locking rings (PEEK), 5 support flanges (PEEK) 2m solvent line (analytical) | 6025.2511 |
| Fuses kit, including: 5 fuses (0.20A, slow-blow, 5 x 20 mm) 15 fuses (overload protection, 2A, slow-blow, 5 x 20 mm) 5 fuses (4A, slow-blow, 6.3 x 32 mm) | 6030.9003 |
| Liquid reservoir for rear seal wash system (2 reservoirs) | 6030.9501 |
| Menu pen | 6300.0100 |
| Maintenance kit for ISO-3100A, LPG-3400A and M, DGP-3600A and M, HPG-3x00 A and M, including: 2 solvent line filters with filter frits, 1 stirrer, ECTFE tubing (1.60 mm OD x 0.75 mm ID), silicone tubing (2.80 mm OD x 1.30 mm ID), PharMed [™] tubing for peristaltic pump, 2 fittings (1/8", PEEK), 2 solvent line locking rings (PEEK), 2 solvent line support flanges (PEEK), 2 knurled head screws (1/4"-28 for ferrule 1/16"), 1 tube connector for 1.0-2.0 mm tube ID, 2 ferrules (1/16" for ¼"-28), 4 ring seals (DR-8), 8 piston seals, 2 mixing chamber gaskets, 4 O-rings (22x2, silicone), 4 pistons (sapphire), 4 filter frits (stainless steel; porosity: 0.5 μm) for inline filter (LPG-3400M and DGP-3600M), 4 filter frits (stainless steel; porosity: 10 μm) for inline filter (ISO-3100A, LPG-3400A, DGP-3600A, HPG-3x00); 6 filter frits for solvent line filter (stainless steel; porosity: 10 μm), 2 piston support hemispheres, 4 piston support rings, 2 micro flow ring seals, 4 valve cartridges (sapphire) for double check valves | 6035.1961 |

| Description | Part No. |
|---|------------------------|
| Maintenance kit for HPG-3200P, including: 1 stirrer, silicone tubing (2.80 mm OD x 1.30 mm ID), PharMed [™] tubing for peristaltic pump, FEP tubing (4.5 mm OD x 3.0 mm ID), 1 tube connector for 1.0- 2.0 mm tube ID, 2 tube adaptor (3.0 mm ID for 1/4"-28), 4 ring seals (DR-8), 8 piston seals, 2 mixing chamber gaskets, 4 pistons (semipreparative), 4 filter frits (stainless steel, porosity: 10 µm) for inline filter, 2 stainless steel solvent line filters, 4 support rings, 4 valve cartridges (sapphire) for double check valves | 6035.1962 |
| Maintenance kit for LPG-3400(AB and MB) and DGP-3600(AB and MB), including: ECTFE tubing (1.6 mm OD x 0.75 mm ID), silicone tubing (2.80 mm OD x 1.30 mm ID), PharMed TM tubing for peristaltic pump, 2 fittings (1/8", PEEK), 2 solvent line locking rings (PEEK), 2 solvent line support flanges (PEEK), 2 knurled head screws (1/4"-28 for ferrule 1/16"), 1 tube connector for 1.0-2.0 mm tube ID, 2 ferrules (1/16" for 1/4"-28), 4 ring seals (DR-8), 8 piston seals, 4 O rings (22x2, silicone), 4 pistons (sapphire), 4 filter frits (titanium; porosity: 2 μ m) for inline filter; 4 filter frits (titanium; porosity: 10 μ m) for inline filter; 2 filter holders (bottom and top parts) for solvent line filter, 6 filter frits for solvent line filter (titanium; porosity: 10 μ m), 2 piston support hemisphere, 4 support rings, 2 micro flow ring seals, 4 valve cartridges (sapphire) for double check valves, 2 mixing chamber gaskets | 6035.1963 |
| Manual injection valve (analytical/micro; \rightarrow section 11.2, page 167) | 5035.0600 |
| Manual injection valve (semipreparative; \rightarrow section 11.2, page 167) | 5035.0601 |
| Micro Flow kit (\rightarrow section 11.2, page 167) for LPG-3400A and DGP-3600A pumps for LPG-3400AB and DGP-3600AB pumps | 6035.1977 6037.1977 |
| Mixing Chamber Extension kit for a total mixing chamber volume of 928 μL (for LPG-3400A, DGP-3600A, HPG-3200A, HPG-3400A pumps) | 6035.1978 |
| Mixing Chamber Extension kit for a total mixing chamber volume of 928 μ L (for LPG-3400AB and DGP-3600AB pumps) | 6037.1978 |
| Mixing Chamber Extension kit for a total mixing chamber volume of 1528 μL (for LPG-3400A, DGP-3600A, HPG-3200A, HPG-3400A, HPG-3200P pumps) | 6035.1979 |
| Mixing Chamber Extension kit for a total mixing chamber volume of 1528 μ L (for LPG-3400AB and DGP-3600AB pumps) | 6037.1979 |
| O-ring seal (22x2; 5 seals for rear seal wash body; all pumps except HPG-3200P) | 6266.2202 |
| Piston (all pumps except HPG-3200P) | 6035.2240 |
| Piston (HPG-3200P) | 2267.0802A |
| Piston support hemisphere (all pumps except HPG-3200P) | 6269.0040 |
| Piston bushing for HPG-3200P (5 bushings) | 2090.8083 |
| Piston retainer for HPG-3200P | 6004.8104 |
| Piston seal/support ring kit (reversed phase, biocompatible) for LPG-3400AB and MB and DGP-3600AB and MB The kit includes 1 support ring and 2 piston seals. | 6025.2012 |
| Piston seal/support ring kit (reversed phase) for ISO-3100A, LPG-3400A or M, DGP-3600A or M, and HPG-3x00A or M. The kit inludes 1 support ring and 2 piston seals. | 6025.2010A |

| Description | Part No. |
|---|------------|
| Piston seal/support ring kit (normal phase) for ISO-3100A, LPG-3400A or M, DGP-3600A or M, and HPG-3x00A or M. The kit includes 1 support ring and 2 piston seals. | 6025.2011A |
| Piston seal/support ring kit (reversed phase) for HPG-3200P. The kit inludes 1 support ring and 2 piston seals. | 6030.9010 |
| Piston seal/support ring kit (normal phase) for HPG-3200P. The kit inludes 1 support ring and 2 piston seals. | 6030.9011 |
| Purge valve | 6020.2030 |
| Rear seal wash body, extractor for | 6000.0046 |
| Rear seal wash system (LPG-3400M and MB, DGP-3600M and MB), including: 1 piston seal), 1 O-ring, 1 rear seal wash body, 1 retainer | 6035.2010 |
| Rear seal wash system (HPG-3200P), including: 1 piston seal, 1 O-ring, 1 rear seal wash body, 1 retainer | 6025.2210A |
| Rear seal wash system (ISO-3100A, LPG-3400A(B), DGP-3600A(B), HPG-3x00A), including: | 6035.2020 |
| Rear seal wash system, tubing kit, including: 1 tube connector 1m silicone tubing Note: Do not use this silicone tubing in the peristaltic pump. For the peristaltic pump tubing, order part no. 6000.5000. | 6030.9502 |
| Retaining guide (solvent bottle) | 6000.0042 |
| Ring seal (DR-8, 10 seals) | 2266.0082 |
| Solvent line (OD 3.0 x ID 1.5 mm) with fittings | 6030.2542 |
| Solvent line from degasser to proportioning valve (DGP-3600A, AB, M, and MB) (set of 3 solvent lines plus line labels) | 6030.2547 |
| Solvent lines from degasser to working head (HPG-3200A, HPG-3200M) (Set of 2 solvent lines plus line labels) | 6035.2530 |
| Solvent lines from degasser to solvent selector (HPG-3400A, HPG-3400M) (Set of 4 solvent lines plus line labels) | 6035.2532 |
| Solvent line (HPG-3200P) | 6250.4301 |
| Solvent reservoir (250 mL) with bottle cap | 2270.0026 |
| Static Mixer kit (for HPG-3x00M pumps) | 6035.0002 |
| Static Mixer - replacement cartridge | 6268.5050 |
| Syringe (100 μ L) for use with manual injection valve (analytical or micro; \rightarrow section 11.2, page 167) | 6035.0665 |
| Syringe (5 mL) for use with manual injection valve (semipreparative; \rightarrow section 11.2, page 167) | 6035.0665 |
| Syringe and tubing kit, including: 5 plastic syringes 3m silicone tubing (2.80 mm OD x 1.30 mm ID) | 6000.0010 |

| Description | Part No. |
|--|-----------|
| T-Insert (HPG-3200 A, HPG-3400 A, and HPG-3200P) | 6035.1980 |
| Tool kit for analytical pumps, including 1 off: Double open-end wrench ($1/4 \ge 5/16$), double open-end wrench ($11 \ge 13$) Allen wrench (size 3), slotted screwdriver (for $1/8$ " pistons) | 6007.9302 |
| Tool kit for semipreparative pump, including 1 off: Double open-end wrench (1/4 x 5/16), double open-end wrench (11 x 13) Allen wrench (size 3), slotted screwdriver (for 8mm pistons) installation tool for cap seal | 6007.9303 |
| Tubing for peristaltic pump (PharMed TM tubing) | 6000.5000 |
| Tubing from proportioning valve to working head (LPG-3400M(B), DGP- 3600M(B)), including the appropriate fittings and ferrules | 6035.2514 |
| Tubing from proportioning valve to working head (LPG-3400A(B) and DGP- 3600A(B)) and from solvent selector to working head (HPG-3400M) including the appropriate fittings and ferrules | 6035.2515 |
| Tubing from degasser to proportioning valve (LPG-3400A, AB, M, or MB) (set of 4) | 6035.2540 |
| USB cable, type A to type B, USB 2.0, 1 m | 6035.9035 |
| USB cable, type A to type B, USB 2.0, 5 m | 6911.0002 |
| Valve cartridge (sapphire) | 6035.2300 |
| Valve nut kit for ISO-3100A, LPG-3400A or M, DGP-3600A or M, and HPG-3x00A or M, including inlet check valve nut and outlet check valve nut | 6035.1965 |
| Valve nut kit for double check valve for LPG-3400AB and MB and DGP-3600AB and MB, including inlet check valve nut and outlet check valve nut | 6037.1964 |
| Valve nut kit for double check valve for HPG-3200P, including: inlet check valve nut and outlet check valve nut | 6035.1966 |

12 Reference Information

12.1 Chemical Resistance of PEEK

For information about the chemical resistance of PEEK, see the table.

- ▲ Important: While PEEK has superb chemical resistance to most organic solvents, it tends to swell when in contact with chloroform (CHCl₃), dimethyl sulfoxide (DMSO), or tetrahydrofuran (THF). In addition, it is attacked by concentrated acids, such as, sulfuric acid and nitric acid or a mixture of hexane, ethyl actate, and methanol. (Swelling or attack by concentrated acids is not a problem with brief flushing procedures.)
- ▲ Important: Bien que PEEK présente une excellente résistance chimique à la plupart des solvants organiques, il a tendance à gonfler lorsqu'il est en contact prolongé avec du chloroforme (CHCl₃), du diméthyle sulfoxyde (DMSO) ou du tétrahydrofurane (THF). De plus, il est attaqué par des acides concentrés tels que l'acide sulfurique et l'acide nitrique ou d'un composé du hexane, éthyle acétate et méthanol. (Ces acides peuvent cependant être utilisés dans le cadre de procédures de nettoyage, à condition que l'exposition soit brève.

| Medium | Concentration [%] | Temperature | Max. Duration | Resistance (+ = yes, - = no) |
|-------------------------|----------------------|-------------|------------------|--|
| Acetaldehyde | techn. pure | 23 | | + |
| Acetone | 100 | 23 | 7 days | + |
| Ammonia | 28 | 23 | 7 days | + |
| Ammonium sulphate | | 23 | | + |
| Amyl acetate | 100 | 23 | | + |
| Amyl alcohol | techn. pure | 23 | | + |
| Benzaldehyde | | 23 | 7 days | + |
| Benzene | 100 | 23 | 7 days | + |
| Benzene/Benzene mixture | | 60 | 42 days | + |
| Benzoic acid | | 23 | | + |
| Borax | | 60 | | + |
| Bromine | | 23 | | - |
| Butane | | 23 | | + |
| Butanol | 100 | 23 | | + |
| Calcium hydroxide | | 23 | | + |
| Carbon dioxide | 100 | 23 | | + |
| Carbon tetrachloride | 100 | 23 | | + |
| Chloric gas | | 23 | | + |

| Medium | Concentration [%] | Temperature | Max. Duration | Resistance (+ = yes, - = no) |
|----------------------|----------------------|-------------|------------------|---------------------------------|
| Chlorine (liquid) | | 23 | | - |
| Chlorobenzene | 100 | 23 | | + |
| Chloroform | 100 | 23 | | + |
| Chromic acid | 40 | 23 | | + |
| Citric acid | | 23 | | + |
| Copper(II) sulphate | | 23 | | + |
| Cyclohexane | 100 | 23 | | + |
| Cyclohexanol | 100 | 23 | | + |
| Cyclohexanone | | 23 | | + |
| Diethyl ether | 100 | 23 | 7 days | + |
| Diisopropyl ether | 100 | 23 | | + |
| Dimethylformamide | 100 | 23 | 7 days | + |
| Dioctylphthalate | | 23 | | + |
| Dioxan | | 23 | | + |
| Ethanoic acid | 96 | 23 | 7 days | + |
| Ethanol | 96 (Vol.) | 23 | 7 days | + |
| Ethyl acetate | 100 | 23 | | + |
| Ethylene glycol | | 23 | | + |
| Ferric chloride | | 23 | | + |
| Formaldehyde | 30 | 23 | | + |
| Formic acid | 95 | 104 | 42 days | + |
| Glycerin | | 23 | | + |
| Heptane | 100 | 23 | 7 days | + |
| Hydrochloric acid | 37 | 23 | | + |
| Hydrofluoric acid | | 23 | | - |
| Hydrogen peroxide | 30 | 23 | 7 days | + |
| Hydrogen sulphide | | 23 | | + |
| Kerosene (Paraffin) | | 23 | | + |
| Lactic acid | | 23 | | + |
| Magnesium chloride | | 23 | | + |
| Methanol | 100 | 23 | | + |
| Methyl ethyl ketone | 100 | 23 | | + |
| Methylisobutylcetone | 100 | 23 | | + |
| Nitric acid | 40 | 23 | 7 days | + |
| Nitric acid | 65 | 23 | 7 days | + |
| Nitrobenzene | 100 | 23 | | + |

| Medium | Concentration [%] | Temperature | Max. Duration | Resistance (+ = yes, - = no) |
|----------------------------|----------------------|-------------|------------------|---------------------------------|
| Paraffin oil | | 60 | | + |
| Perchlorethylene | 100 | 23 | | + |
| Phenol | diluted | 23 | | + |
| Phenol | conc. | 23 | | - |
| Potassium dichromate | | 23 | | + |
| Potassium hydroxide | | 23 | | + |
| Potassium nitrate | | 23 | | + |
| Potassium permanganate | | 23 | | + |
| Propane | | 23 | | + |
| Propyl alcohol | | 100 | | + |
| Pure benzene | | 60 | | + |
| Silicone oil | | 160 | | + |
| Sodium chloride | | 23 | | + |
| Sodium hydrogen carbonate | | 23 | | + |
| Sodium hydroxide | 40 | 23 | 7 days | + |
| Sodium hydroxide | 30 | 130 | | + |
| Sodium thiosulphate | | 23 | | + |
| Sulphur dioxide | | 23 | | + |
| Sulphuric acid | 40 | 130 | | + |
| Sulphuric acid | 50 | 23 | 7 days | + |
| Sulphuric acid (dissolved) | 98 | 23 | | - |
| Toluol | 100 | 23 | 7 days | + |
| Trichloroethylene | 100 | 23 | 7 days | + |
| Water | | 23 | | + |
| Xylene | 100 | 23 | | + |
| Zinc chloride | | 23 | | + |

12.2 Solvent Miscibility

Miscibility describes the ability of liquids to form homogeneous mixtures in all proportions (one-phase system). Solvent miscibility is important during elution and when changing from one solvent to another. Thus, when you prepare solvents consider the miscibility and homogeneous mixing of the single components. Note that certain compositions of some solvent systems may result in miscibility gaps.

For information about solvent miscibility, see the table (source: *Handbuch der HPLC*, GIT Verlag, 1995). The table provides a general idea of solvent miscibility. Under certain conditions, non-miscible liquids may mix or miscible liquids may separate.



12.3 Properties of Common Solvents

The table summarizes the properties of the most important solvents in HPLC [1, 2]

| | Acetronitrile | Dichloromethane | n-Hexane | Isorpopanol | Methanol | Tetrahydrofurane | Water |
|--|---------------|-----------------|----------|-------------|----------|------------------|-------|
| UV Transmission at [nm] | | | | | | | |
| 20% (0.7 AU) | 190 | 235 | 200 | 210 | 210 | 255 | |
| 80% (0.1 AU) | 195 | 245 | 225 | 230 | 235 | 370 | |
| 98% (0.01 AU) | 220 | 260 | 260 | 260 | 260 | 310 | < 190 |
| Refraction Index (RI) at 20 °C | 1.344 | 1.424 | 1.376 | 1.378 | 1.329 | 1.406 | 1.333 |
| Boiling Point (BP) in °C at 1013 hPa | 82 | 40 | 69 | 82 | 65 | 66 | 100 |
| Vapor Pressure (VP) at 25 °C | 118 | 582 | 202 | 60 | 169 | 216 | 32 |
| Viscosity (η) at 20 °C (cP = mPa*s) | 0.37 | 0.44 | 0.33 | 2.3 | 0.60 | 0.55 | 1.00 |
| Density (p) (g/mL) | 0.78 | 1.32 | 0.66 | 0.78 | 0.79 | 0.88 | 0.997 |
| $\eta/\rho (cP*mL/g)$ | 0.47 | 0.33 | 0.50 | 2.9 | 0.76 | 0.62 | 1.00 |
| Compressibility (χ) at 20 °C (Mbar ⁻¹) | 99 | 97 | 160 | 100 | 123 | 93 | 46 |
| Critical Flow F _c (mL/min) ¹⁾ | 13 | 9.4 | 14 | 83 | 21 | 18 | 28 |
| Linear Drop in Pressure $\Delta p/l$ (MPa/m) ²⁾ | 0.06 | 0.08 | 0.06 | 0.40 | 0.10 | 0.10 | 0.17 |
| Polarity (P') ³⁾ | 5.8 | 3.1 | 0.1 | 3.9 | 5.1 | 4.0 | 10.2 |

¹⁾ F_C = critical flow for 0.25 mm ID tubing

 F_{C} (mL/min) = 113 x 0.25 mm x η (cP) / ρ (g/mL)

Fc is an example of a hydrodynamic calculation.

²⁾ $\Delta p/l =$ linear drop in pressure for 1 mL/min and 0.25 mm ID tubing $\Delta p/l (MPa/m) = 6.8 \times 10^{-6} \times 1 \text{ mL/min} \times 100 \text{ cm} \times \eta (cP) / (0.25 \text{ mm})^4$ $\Delta p/l$ is an example of a hydrodynamic calculation.

³⁾ P' is the polarity calculated by L.R. Snyder [3] from experimental measurements by L. Rohrschneider [4].

References

- [1] K.K. Unger, E. Weber (Hrsg.), Handbuch der HPLC, GIT Verlag, 1995
- [2] D.R. Lide, Handbook of Chemistry and Physics, 79th Edition, CRC Press, 1998-1999
- [3] L.R. Snyder, Journal of Chromatographic Sciences, 16, 223, 1978
- [4] L. Rohrschneider, Analytical Chemistry, 45, 1241, 1973

12.4 Safety Information about Flammable Solvents

The following table provides an overview of safety information for flammable solvents in HPLC

| | Acetonitrile | Diethylether | Ethanol | Ethylacetate | Heptane | Hexane | Isopropanol | Methanol | Tetrahydofurane |
|--------------------------------|--------------|--------------|---------|--------------|---------|--------|-------------|----------|-----------------|
| Boiling point (°C) | 82 | 35 | 78 | 77 | 98 | 69 | 82 | 65 | 66 |
| Vapor pressure (hPa) | 118 | 735 | 93 | 121 | 55 | 202 | 60 | 169 | 216 |
| Flash point (°C) | 6 | -45 | 12 | -4 | -4 | -22 | 12 | 11 | -14 |
| Auto-ignition temperature (°C) | 520 | 190 | 490 | 490 | 230 | 260 | 540 | 510 | 320 |
| Explosion Limits (%) | 3-16 | 2-36 | 3-19 | 2-36 | 1-7 | 1-8 | 2-12 | 7-36 | 2-12 |

The table is based on the following definitions and references.

Definitions

- The flash point is the lowest temperature at an atmospheric pressure of 1013 mbar at which a liquid gives off enough vapors to ignite with an external ignition source when mixing with the air above the liquid. [1]
- Substances whose flash point is below 38 °C are classified as flammable. [2]
- The auto-ignition temperature is the lowest temperature at which substances can selfignite at atmospheric pressure without an external ignition source, that is, without external ignition by sparks or flames. The thermal energy required to reach the autoignition temperature is created by a spontaneous chemical reaction or physical processes in or on the surface of the combustible substances. The determination of the auto-ignition temperature is imprecise and depends on the equipment and apparatus in use. Nevertheless, it indicates the maximum permissible surface temperature of equipment and apparatus when they are exposed to an air-vapor mixture of these substances. [1, 2]
- The explosion limit are the upper and lower concentration limits of a mixture of a flammable gas or vapor with air in which this mixture can explode when being heated or by means of sparks. [1]

i Tip: Volatile solvents are not necessarily flammable as well. For example, chloroform is volatile but non-flammable.

References

- [1] Otto-Albrecht Neumüller, Römpps Chemie-Lexikon, 8. Auflage, 1987
- [2] W.E. Baker et al., Explosion Hazards and Evaluation, Elsevier Sci. Publ., 1983
- [3] H. Bennett, Concise Chemical and Technical Dictionary, Edward Arnold Ed., 1986
- [4] D.R. Lide, Handbook of Chemistry and Physics, 79th Edition, CRC Press, 1998-1999
- [5] G.W.C. Kaye and T.H. Laby, *Tables of Physical and Chemical Constants*, 16th Edition, Longman Ed., 1995
- [6] Union des Industries Chimiques, *L'Electricité Statique en Atmosphère Explosive*, Septembre 1982
- [7] B.P. Mullins, Spontaneous Ignition of Liquid Fuels, Butterworths Ed., 1955
- [8] Chemical Safety Sheets, Working Safely with Hazardous Chemicals, Kluwer Acad. Publ., Samson Chem. Publ., Dutch Inst. for the Working Environment, and Dutch Chem. Ind. Assoc., 1991
- [9] F.A. Williams, Combustions Theory, Benjamin / Cummings Publ., 1985 Technischer Anhang

13 Appendix

13.1 Connections in the Low- and High-Pressure Sections

The illustration shows the solvent connections in the low-pressure section of the pump.



| No. | Description |
|-----|-------------|
| 1 | Ferrule |
| 2 | Fitting |
| 3 | Capillary |

Fig. 75: Solvent line connection on low-pressure side

Do not overtighten these fitting connections; the connections should be no more than handtight. Tighten more if a connection is leaking. Avoid cross-threading when installing the fittings on the PTFE valve block. Cross-threading might damage the valve blocks.

All capillary connections in the high-pressure section of the pump include fitting screws and ferrules.



Fig. 76: Fitting screw and ferrule

Do not overtighten these fitting connections; the connections should be hand-tight plus an additional one-quarter turn. Tighten more if a connection is leaking. To connect the capillaries to an injection valve or selector valve, install only the ferrules and fittings shipped with the valve and observe the manufacturer's installation instructions.

13.2 Manual Injection Valve

A manual injection value is available as an option for the pump. The mounting angle provides also a storage location for the menu pen.



Fig. 77: Manual injection valve mounted to the pump enclosure

| Description | Part No. |
|---|-----------|
| Manual injection valve (analytical/micro) | 5035.0600 |
| Manual injection valve (semipreparative) | 5035.0601 |

I Tip: The following syringes are available for use with the manual injection valve:

| Description | Part No. |
|---|-----------|
| 100 µL syringe for analytical/micro injection valve | 6035.0665 |
| 5 mL syringe for semipreparative injection valve | 6035.0670 |

If you use a syringe other than those mentioned above, only use syringes with the correct needle (outer diameter 0.028", length 2").

Attach the valve to the bottom right of pump enclosure as described in the installation instructions shipped with the valve. Set up the valve in Chromeleon.

- 1. Start the Chromeleon Server Monitor if it is not already running (\rightarrow page 34).
- 2. Start the Chromeleon Server Configuration program by clicking Start on the taskbar. Point to Programs (or All Programs, depending on the operating system), point to Chromeleon, and then click Server Configuration.
- 3. Select the timebase to which the injection valve will be assigned.

- 4. Open the Add device to timebase dialog box. To do so, click Add Device on the Edit menu or right-click the timebase and click Add Device on the menu.
- 5. On the **Manufacturers** list, click **Generic** and on the **Devices** list, click **Remote Inject**.
- 6. On the **General** page, select a digital input of the pump on the **Inject Port** list, for example, select UM3PUMP_Input1.

| Remote Inject | | × |
|---|-------------|---|
| General BCD Position Inp | outs | |
| | | 1 |
| Device <u>N</u> ame: | InjectValve | |
| Inject Port: | | |
| | | |
| <none2< td=""><td></td><td></td></none2<> | | |
| UM3PUMP Input1 | | |
| UM3PUMP_Input2 | | |
| UM3PUMP_Input3 | | |
| TCC3x00_INPUT_1 | | |
| [TCC3x00_INPOT_2 | | |
| | | |
| | | |
| | | |
| | | |
| | | |

Fig. 78: Setting up the valve for external injection

The manual injection valve now available in Chromeleon and appears in the Chromeleon **Commands** dialog box under **InjectValve**.

- 7. Add an **Inject** command to the Chromeleon program. If you create the program with the Program Wizard, the **Inject** command is added automatically.
- 8. When Chromeleon executes the program, it runs the commands that occur before the **Inject** command, and then waits for the signal from the injection valve. The following message is displayed in the Audit Trail "Wait for inject response on remote start." When the inject signal occurs, Chromeleon resumes execution of the program.

13.3 Pin Assignment

13.3.1 D-Sub I/O (Digital I/O)

| Pin | Wire | Signal Name | Signal Level | Remarks |
|-----|------|-------------|----------------|------------------------------|
| 1 | 1 | | | Marked wire/reserved |
| 2 | 3 | | | Reserved |
| 3 | 5 | RELAY 3 OUT | Potential free | Closing contact |
| 4 | 7 | RELAY 1 OUT | Potential free | Opening contact |
| 5 | 9 | RELAY 2 OUT | Potential free | Opening contact |
| 6 | 11 | RELAY 3 OUT | Potential free | Opening contact/Operable Out |
| 7 | 13 | RELAY 1 OUT | Potential free | Middle contact |
| 8 | 15 | RELAY 2 OUT | Potential free | Middle contact |
| 9 | 17 | GND | Ground | Reference potential |
| 10 | 19 | HOLD IN | Ground | Reference potential |
| 11 | 21 | STOP IN | Ground | Reference potential |
| 12 | 23 | START IN | Ground | Reference potential |
| 13 | 25 | | | Reserved |
| 14 | 2 | RELAY 4 OUT | Potential free | Closing contact |
| 15 | 4 | RELAY 4 OUT | Potential free | Middle contact |
| 16 | 6 | RELAY 4 OUT | Potential free | Opening contact |
| 17 | 8 | | | Reserved |
| 18 | 10 | RELAY 3 OUT | Potential free | Middle contact/Operable Out |
| 19 | 12 | RELAY 1 OUT | Potential free | Closing contact |
| 20 | 14 | RELAY 2 OUT | Potential free | Closing contact |
| 21 | 16 | Vcc_Save | +5V/500mA | |
| 22 | 18 | HOLD IN | TTL | Digital input 1 |
| 23 | 20 | STOP IN | TTL | Digital input 2 |
| 24 | 22 | START IN | TTL | Digital input 3 |
| 25 | 24 | | | Reserved |

Fig. 79: 25-pin D-Sub I/O port (female)

 \triangle Important: The maximum switching voltage of the relays is 24 V. The switching current must not exceed 100 mA.

Important: La tension maximale de commutation des relais est de 24 V. L'intensité de commutation ne doit pas dépasser 100 mA.

| Pin | Signal Name | Signal Level | Remarks |
|-----|--------------------|---------------|-----------------------------------|
| 1 | | | Reserved |
| 2 | Solvent Rack Error | | TTL_high with solvent rack errors |
| 3 | | | Jumper to pin 9 |
| 4 | Solvent Rack Leak | | TTL high with solvent rack leaks |
| 5 | | | Reserved |
| 6 | V_Degas | +24V_supply | Supply for the solvent rack |
| 7 | GND_Degas | Ground_supply | Reference potential for VC_Degas |
| 8 | VCC | +5V | Voltage for logic devices |
| 9 | | | Jumper to pin 3 |
| 10 | GND | | Reference potential for VCC |
| 11 | GND | | Reference potential for VCC |
| 12 | GND | | Reference potential for VCC |
| 13 | | | Reserved |
| 14 | V_Degas | +24V_supply | Supply for the solvent rack |
| 15 | GND_Degas | Ground_supply | Reference potential for VC_Degas |

13.3.2 Solvent Rack

Fig. 80: 15-pol. Solvent Rack port (female)

14 Index

A

| Accessories | |
|--------------------------|--|
| Optional | |
| Standard | |
| Active Rear Seal Washing | |
| Actual Operating Hours | |

B

| Back | 69 |
|---------------------|----|
| Batch | 66 |
| Biocompatible Pumps | 14 |
| Bootloader version | 73 |

С

| | 69 |
|-------------------------------------|----|
| Check Valve Replacement | 13 |
| Chromeleon | 61 |
| Automated Control | 65 |
| Connect | 61 |
| Diagnostics | 27 |
| Direct Control | 62 |
| Installation Program | 34 |
| License | 26 |
| Load USB Driver | 34 |
| Predictive Performance | 86 |
| Program Wizard | 65 |
| Pump Configuration | 37 |
| Pump Installation | 36 |
| Server Configuration Program | 36 |
| Server Monitor | 36 |
| Special Functions | 86 |
| Chromeleon Dialog Boxes | |
| Configuration complete | 41 |
| Devices | 39 |
| Error Levels | 44 |
| General | 37 |
| Head Type & Limits | 42 |
| Inputs | 43 |
| Pump (Right/Left) | 42 |
| Relays | 43 |
| Solvents | 43 |
| Solvents (Right/Left) | 43 |
| Cleaning (Piston) | 24 |
| Column Flow | 60 |
| Commands (Dialog Box) | 62 |
| Commands (Display) | 67 |
| Compression | 73 |
| Configuration complete (Dialog Box) | 41 |
| Configuration Menu | 74 |
| Configurations | 12 |

| Connection | |
|-----------------------|--|
| Chromeleon Computer | |
| Power Cord | |
| USB | |
| Connections | |
| Fluid | |
| High-Pressure Section | |
| Low-Pressure Section | |
| Consumables | |
| Control (Automated) | |
| Control (Direct) | |
| Control Elements | |
| Control Menu | |
| Control Panel | |
| | |

D

| DAC Output | 74 |
|----------------------|-------------|
| Degasser | 72 |
| Device View | 66 |
| Devices (Dialog Box) | 39 |
| DGP-3600 | 148 |
| Fluid Connections | 149 |
| Inside Front Panel | 148 |
| Diagnostics | |
| Diagnostics | |
| Diagnostics Menu | 73 |
| Diagnostics Messages | |
| Digital I/O | 20, 88, 186 |
| Display | |
| Display & soft keys | 74 |
| Display Mode | 74 |
| Display Mode) | 77 |
| Double Flow Mode | 43, 156 |
| Double Flow Mode) | 89 |
| Drain System | 53 |
| Drainage | |
| Pump | |
| System | |
| | |

Е

| Equilibration | 57 |
|-------------------------------------|----|
| Equilibration (SmartStartup Wizard) | 57 |
| Error Levels (Dialog Box) | 44 |
| Error Messages | |

F

| Facility Requirements | |
|-----------------------------|------------|
| Filter Frit (Inline Filter) | |
| Firmware Download | |
| Firmware Version | 73 |
| Flow Acceleration | 72, 78, 79 |

| Flow Deceleration | .72,78 | 8, 79 |
|---------------------------|--------|-------|
| Flow Off | 60 | 0, 67 |
| Flow On | 6 | 0, 67 |
| Flow Rate | | 78 |
| Flow Splitter | | 77 |
| Fluid Connections | | 21 |
| DGP-3600 | | . 149 |
| High-Pressure Section | | .183 |
| HPG-3200 | | .151 |
| HPG-3400 | | .154 |
| ISO-3100 | | .145 |
| Low-Pressure Section | | .183 |
| LPG-3400 | | .147 |
| Rear Seal Washing | | 49 |
| Fluid Connections | | .143 |
| Found New Hardware Wizard | | 35 |
| Front Panel | | 17 |
| Front Panel Display | | |
| Function Key | | 67 |
| Power-Up | | 59 |
| Screen Brightness | | 85 |
| Screen Contrast | | 85 |
| Status Screen | | 60 |
| Function Key | | 67 |
| Arrow down | | 69 |
| Arrow right | | 69 |
| Arrow up | | 69 |
| Back | | 69 |
| Cancel | | 69 |
| Clear | 7(| 0, 96 |
| Flow Off | | 67 |
| Flow On | | 67 |
| Menu | | 67 |
| Next | 7(| 0, 96 |
| Ok | | 69 |
| Prev | 70 | 0, 96 |
| Purge | | 67 |
| Select | | 69 |
| Set Flow | | 67 |
| Fuse Cartridge | | 19 |
| Fuse Replacement | | .131 |

G

| General (Dialog Box) | |
|-----------------------|-----|
| Gradient Combinations | 156 |
| Gradient Curve | |

H

| Head Type & Limits (Dialog Box) | |
|-------------------------------------|-----|
| High-Pressure Gradient | 156 |
| Double Flow Mode | 156 |
| Gradient Combinations | 156 |
| Semipreparative Pump | 157 |
| High-Pressure Gradient Pump | 150 |
| High-Pressure Section (Connections) | |

| HPG-3200 | |
|--------------------|-----|
| Fluid Connections | 151 |
| Inside Front Panel | 150 |
| HPG-3400 | |
| Fluid Connections | |
| Inside Front Panel | |

I

| Inject Synchronization | |
|------------------------|---------|
| Inline Filter | 23, 111 |
| Exchange | 112 |
| Filter Frit | 112 |
| Inputs (Dialog Box) | |
| Inside Front Panel. | 143 |
| DGP-3600 | 148 |
| HPG-3200 | 150 |
| HPG-3400 | 153 |
| ISO-3100 | 144 |
| LPG-3400 | 146 |
| Installation | |
| Connection | |
| Digital I/O | |
| in Chromeleon | |
| Location | |
| Power Connection | |
| Sample Injector | 184 |
| Solvent Rack | |
| Instrument Shut Down | |
| Intended Use | |
| Interior Components | |
| ISO-3100 | |
| Fluid Connections | 145 |
| Inside Front Panel | 144 |
| Isocratic Pump | 144 |

L

| Leak Detection | |
|------------------------------------|-------------|
| Leak Rate | |
| Leak Sensor | 25, 84, 110 |
| Leak Sensor Mode | |
| Leakage Test | |
| Left Pump (Dialog Box) | |
| Left Solvents (Dialog Box) | |
| Linking | |
| Autosampler-Pump | |
| Pump-Flow Manager | 77 |
| Low-Pressure Gradient Pump | |
| Low-Pressure Section (Connections) | |
| LPG-3400 | |
| Fluid Connections | |
| Inside Front Panel | |
| Vacuum Degasser | 25, 132 |

Μ

| Main Menu | |
|--------------------------|-----|
| Maintenance | |
| Maintenance Intervals | |
| Master Flow | 60 |
| MaximumFlowRampDown | |
| MaximumFlowRampUp | |
| Menu | |
| Configuration | 74 |
| Control | 71 |
| Diagnostics | 73 |
| Layout | 69 |
| Main | |
| Overview | |
| Preferences | 72 |
| Structure | 69 |
| Messages | 96 |
| Micro Flow Kit | |
| Mixer (Static) | 141 |
| Mixing Chamber | |
| Mixing Chamber Extension | |
| Model | 73 |
| Module Address | |
| | |

0

| Off | 60 |
|---------------------------|---------|
| Ok | 69 |
| On | 60 |
| Operating Instructions | 1 |
| Operation | 59 |
| Chromeleon | 61 |
| Control Elements | 17 |
| Front Panel Display | 17 |
| Front Panel Keys | 67 |
| Function Key | 67 |
| Power-Up | |
| Safety Precautions | 3 |
| Safety Symbols | 2 |
| Solvent Selection | 75, 175 |
| Operation Principle | |
| Operational Qualification | |
| Operational Settings | 75 |
| Optimizing | 135 |
| Micro Flow Kit | 138 |
| Mixing Chamber Extension | 136 |
| Static Mixer | 141 |
| Outlet Unit | 11.112 |
| Inline Filter | |
| Mixing Chamber | 24 |

Р

| Panel Tabset | 64 |
|---------------------------|-----|
| Performance Qualification | 89 |
| Pin Assignment | 186 |

| Piston | |
|------------------------|--------|
| Install | 125 |
| Remove | 120 |
| Piston Cleaning | 124 |
| Piston Seal Washing | 22, 49 |
| Alternative Setup | |
| Piston Seals | 114 |
| Inspecting for Leakage | |
| Install | 125 |
| Remove | 123 |
| Replace | 119 |
| Power Connection | |
| Power Switch | 19 |
| Power-Up | |
| Predictive Performance | 27, 86 |
| Preferences Menu | |
| Preparation | |
| Actions | |
| Capillaries | |
| Equilibration | |
| Pressure Limits | 79 |
| Pressure Output | |
| Pressure Unit | 74 |
| Principle of Operation | 10 |
| Problems | 103 |
| Program Start | |
| Program Wizard | 65 |
| Pump (Dialog Box) | |
| Pump Design | 16 |
| Pump Diagnostics | |
| Pump Drainage | 53 |
| Pump Head | |
| Install | 125 |
| Remove | 120 |
| Pump Pressure | 80 |
| Purge | 60, 67 |
| Purge Flow | 72 |
| Purge Time | 72 |
| Purging | 54, 84 |
| Manually | |
| Standard Settings | 54 |
| Via the Autosampler | |

R

| Rear Panel | |
|---------------------------|------------|
| Rear Panel Connector | |
| Digital I/O | |
| Pressure Output | |
| Solvent Rack | |
| Rear Seal Limit | 72 |
| Rear Seal Wash | |
| Rear Seal Washing | 22, 49, 72 |
| Alternative Setup | |
| Connect | |
| External Liquid Reservoir | |
| | |

| Internal Liquid Reservoir | 50 |
|-----------------------------|-----|
| Standard Setup | 22 |
| Turn on | 81 |
| Washing Medium | 51 |
| Relays (Dialog Box) | 43 |
| Reset to Factory Defaults | 74 |
| Right Pump (Dialog Box) | 42 |
| Right Solvents (Dialog Box) | 43 |
| RMA Number | 109 |

S

| Safety Information | 2 |
|-------------------------------------|---------|
| Safety Precautions | 3 |
| Safety Symbols | 2 |
| Sample Injector (Installation) | |
| Screen Brightness | |
| Screen Contrast | |
| Seal Washing | 22, 49 |
| Select | 69 |
| Selection (Solvent) | 75, 175 |
| Self Test. | 73 |
| Semipreparative Pump | 157 |
| Serial number | 73 |
| Server Configuration Program | |
| Server Monitor | |
| Service | |
| Check Valve Replacement | |
| Filter Frit (Inline Filter) | |
| Filter Frit (Outlet Unit) | |
| Fuse Replacement | |
| General Notes | |
| Inspecting Piston Seals for Leakage | 117 |
| Leak Sensor | |
| Vakuum Degasser | |
| Settings (Operation) | 75 |
| Shut Down | 90 |
| Shutdown Program | 91 |
| SmartShutdown | 85, 91 |
| SmartStartup | 57, 85 |
| Soft Key | 67 |
| Software Control | 26, 61 |
| Solvent Lines | |
| Solvent Rack | 21, 187 |
| Solvent Reservoirs | 47 |
| Connect | |
| General Precautions | 47 |
| Solvents | |
| Chemical Resistance (PEEK) | |
| Miscibility | |
| | |

| Properties | |
|--------------------------|--|
| Safety Information | |
| Selection | |
| Solvents (Dialog Box) | |
| Spare Parts | |
| Specification | |
| Standby Program | |
| Startup | |
| Actions | |
| Capillaries | |
| Equilibration | |
| Solvent Lines | |
| Static Mixer Kit | |
| Status Screen | |
| Sync Inject | |
| Synchronization (Inject) | |
| System Wellness | |

Т

| Technical Information | |
|-----------------------|----------|
| Tightness | 110, 117 |
| Total Operating Hours | |
| Troubleshooting | |
| Diagnostics Messages | |
| Messages | |
| Operating Problems | |

U

| Unit Description | 9 |
|------------------------|---|
| Biocompatible Pumps | |
| Configurations | |
| Overview | 9 |
| Principle of Operation | |
| Semipreparative Pump | |
| Unpacking | |
| USB | |
| User Information | |
| | |

V

W

| Wash Liquid (Degas) | 4′ | 7 |
|---------------------|----|---|
| Workload | 7. | 3 |