

Thermo Scientific Dionex UltiMate 3000 Series

SR-3000 Solvent Rack

Operating Instructions (Original Operating Instructions)



Revision: 1.7

Date: September, 2013

© 2013 Thermo Fisher Scientific Inc.
Doc. No. 4820.3555





Declaration of Conformity

(Original Declaration of Conformity)

Product: Thermo Scientific Dionex UltiMate 3000 Solvent Rack

Types: **SR-3000, SRD-3200, SRD-3400, SRD-3600**

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

- Low-Voltage Directive 2006/95/EC
- EMC Directive 2004/108/EC

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

- DIN EN 61010-1:2010
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

This declaration is issued for the manufacturer

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

by the Managing Director, Rüdiger Obst and
the Vice President HPLC, Fraser McLeod.

September 2, 2013

Table of Contents

1	Introduction.....	1
1.1	How to Use This Manual	1
1.2	Safety Information	2
1.2.1	Symbols on the Instrument and in the Manual	2
1.2.2	Safety Precautions.....	3
1.2.3	Consignes de Sécurité.....	6
1.3	Intended Use	9
1.4	Federal Communications Commission (FCC) Note.....	9
2	Installation and Maintenance	11
2.1	General Information.....	11
2.2	Unpacking	11
2.3	Positioning the Solvent Rack in the UltiMate 3000 System.....	12
2.4	Inside View	13
2.5	Solvent Reservoirs	14
2.5.1	General Precautions	14
2.5.2	Connecting the Solvent Reservoirs.....	15
2.6	Routine and Preventive Maintenance	16
3	Technical Information.....	17
4	Accessories, Consumables, and Spare Parts	19
4.1	Standard Accessories	19
4.2	Consumables and Spare Parts	19

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 Thermo Scientific Dionex™ UltiMate™ 3000 Solvent Rack, Thermo Fisher Scientific recommends that you review the manual thoroughly before beginning operation.

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

PEEK is a trademark of Victrex PLC.




All other trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries.

1.2 Safety Information






The CE Mark label on the instrument indicates that the instrument is compliant with the related standards.

1.2.1 Symbols on the Instrument and in the Manual

The table shows the symbols used on the instrument:


Symbol	Description
	WEEE (Waste Electrical and Electronic Equipment) label—For more information, see the WEEE Information section in the "Installation and Qualification Documents for Dionex Thermo Scientific Instruments" binder. Étiquette DEEE (Déchets d'Equipements Electriques et Electroniques) — Pour plus d'informations, référez-vous au chapitre WEEE Information dans le classeur "Installation and Qualification Documents for Chromatography Instruments".
	Refer to the operating instructions 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.
	Label according to the "Measures for Administration of the Pollution Control of Electronic Information Products" (China RoHS) guideline if necessary Étiquette China RoHS si nécessaire

At various points throughout the manual, the following symbols indicate messages of particular importance:


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

1.2.2 Safety Precautions

When working with analytical instrumentation, you must know the potential hazards of using chemical solvents.

 **Tip:** Before initial operation, make sure that you are familiar with the contents of this manual.

For the safety precautions in French, see section 1.2.3 (→ page 6).

 **Warning:** All users of the device must observe the following safety precautions and all additional safety precautions in this manual to avoid the possibility of personal injury or damage to the device when operating the device or carrying out any maintenance or service procedures.

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

- **Protective equipment**

When performing any work on or near the HPLC system, wear personal protective equipment (protective clothing, safety gloves, safety glasses) as required by the hazard of the mobile phase and sample. For information about the proper handling of particular substance and for advice on specific hazards, refer to the material safety data sheet for the substance you are using. Observe the guidelines of Good Laboratory Practice (GLP).

An eyewash facility and a sink should be close to the device. If any substance splashes on the eyes or skin, wash the affected area and seek medical attention.

- **Hazardous substances**

Many organic solvents, mobile phases, and samples are harmful to health. Be sure that you know the toxic and infectious properties of all substances that you are using. You may not know the toxic or infectious properties of many substances that you are using. If you have any doubt about a substance, treat it as if it contains a potentially harmful substance. For advice on the proper handling of a particular substance, refer to the Safety Data Sheet (SDS) of the manufacturer. Observe the guidelines of Good Laboratory Practice (GLP).

Dispose of waste substance in an environmentally safe manner that is consistent with all local regulations. Do not allow flammable, toxic, and/or infectious substances to accumulate. Follow a regulated, approved waste disposal program. Never dispose of flammable, toxic, and/or infectious substances through the municipal sewage system.

- **Hazardous gases**

Install the HPLC system in a well-ventilated laboratory. If the mobile phase or sample includes volatile or flammable solvents, do not allow them to enter the workspace. If the mobile phase or sample includes volatile or flammable solvents, avoid open flames and sparks.

- **Electrostatic discharge**

Discharge of electrostatic energy may lead to sparking and can constitute a fire hazard. Keep in mind that liquid flowing through capillaries can generate static electricity. This effect is particularly pronounced in insulating capillaries and with non-conductive solvents (for example, pure acetonitrile).

Take appropriate measures to prevent the generation of static electricity near the HPLC system. For example, make sure that the air humidity level in the laboratory is sufficiently high and provide proper ventilation, wear anti-static clothing or shoes, prevent accumulation of air bubbles in waste lines, and use grounded waste containers. Use only non-conductive capillaries to direct solvents into the waste container. With electrically conductive capillaries, make sure that they are properly grounded.

- **Self-ignition of solvents**

Do not use solvents for which the self-ignition temperature is below 150 °C. In case of leakage, these solvents may self-ignite on a hot surface.


- **Capillaries, capillary connections, open connections**


- ◆ Capillaries, especially non-metallic capillaries may burst, slip out of their fittings or may not be screwed in. This may result in substances spraying out of the open connections.
 - ◆ In an UltiMate 3000 system, some components are made of PEEK™. This polymer has superb chemical resistance to most organic solvents. However, it tends to swell when in contact with trichloromethane (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 acetate, and methanol. In both cases, capillaries may start leaking or they can burst. Swelling or attack by concentrated acids is not a problem with brief flushing procedures.
 - ◆ Do not use tubing that is stressed, bent, kinked, or damaged.
 - ◆ Capillary connections can be contaminated by harmful substances or harmful substances can escape from open connections.
 - ◆ Some capillaries of the RS pumps and some Viper system capillaries are made of MP35N[®], a nickel-cobalt based alloy. Individuals with sensitivity to nickel/cobalt may show an allergic reaction from skin contact.
 - ◆ Always wear safety glasses when handling fused silica tubing, for example, during installation or when cutting capillaries to the length.
- For safety reasons, you shall not lift the SRD with one or more solvent reservoirs being in the tray.
 - Use only the original spare parts and accessories authorized for the device by Thermo Fisher Scientific.

- When operating the HPLC system, always set a lower pressure limit for the pump. This prevents damage resulting from leakage or from running the pump dry over a longer period. Activate solvent reservoir level monitoring (→ *Pump manual*).
- The front panel tilts upward. When lifting or moving the SRD, always lift by the bottom or sides of the instrument. This is to avoid damage to the instrument.
- The open front panel door is not designed to carry weight. Therefore, you shall not place any objects on the open front panel door.
- Do not use the solvent rack in ways other than those described in this manual.
- Keep the operating instructions near the device to be available for quick reference.

1.2.3 Consignes de Sécurité

Si vous utilisez d'instrumentation analytique, vous devez connaître les risques d'utilisation de produit chimiques.

 **Veillez noter:** Avant de commencer à utiliser l'instrument, assurez-vous que vous vous êtes familiarisés avec le contenu de ce manuel.

 **Avertissement:** Toutes les personnes utilisant l'instrument doivent observer les consignes de sécurité suivantes et dans les autres chapitres de ce manuel pour éviter une mise en danger de sa personne ou de dommage à l'instrument pendant l'utilisation et des opérations de maintenance ou service de l'instrument.

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

- **Equipment de protection**

Pour tous les travaux sur le système HPLC ou à proximité, portez l'équipement de protection personnel (vêtements de protection, gant de sécurité, lunettes de protection) qui correspond aux risque découlant de la phase mobile et/ou de l'échantillon. Pour les informations sur la manipulation correcte des composés et des recommandations pour les situations de risque spécifiques, veuillez consulter la fiche de données de sécurité des substances que vous utilisez. Veuillez respecter des directives des Bonnes Pratiques de Laboratoire (BPL).

Une installation permettant de se laver les yeux ainsi qu'un lavabo doivent se trouver à proximité du système. Si une substance, quelle qu'elle soit, entre en contact avec vos yeux ou votre peau, rincez abondamment la zone affectée à l'eau, puis.

- **Substances dangereuses**

De nombreux solvants organiques, phases mobiles et échantillons sont nuisibles à la santé. Informez-vous de propriétés toxicologiques et infectieuses de toutes les substances que vous utilisez. Les propriétés toxicologiques et infectieuses de nombreuses substances peuvent être mal connues. Au moindre doute concernant une substance, traitez-la comme s'il contenait une substance potentiellement dangereuse. Pour des instructions comment utiliser correctement des composés particuliers, veuillez consulter à la fiche de données des sécurités du fabricant respectif. Veuillez respecter des directives des Bonnes Pratiques de Laboratoire (BPL).

Débarressez-vous de tous les déchets de substances de manière écologique, conformément à la réglementation en vigueur au niveau local. Empêchez impérativement l'accumulation de solvants inflammables, toxiques et/ou infectieux. Suivez un programme d'élimination des déchets réglementé et approuvé. Ne jetez jamais de solvants inflammables, toxiques et/ou infectieux dans le système municipal d'évacuation des eaux usées.

- **Gaz dangereux**

Installez le système HPLC dans un laboratoire bien ventilé. Si la phase mobile ou l'échantillon contient des solvants volatils ou inflammables, vous devez assurer qu'ils ne pénètrent dans l'espace de travail. Si la phase mobile ou l'échantillon contient des solvants volatils ou inflammables, évitez les flammes nues et les sources d'étincelles à proximité.

- **Décharge électrostatique**

Décharge électrostatique peut provoquer la formation d'étincelles et peut présenter un risque d'incendie. Veuillez noter que des solvants fluides dans les capillaires peuvent se charger automatiquement. Cet effet se peut produire particulièrement forte dans les capillaires isolants et avec des solvants non-conducteurs (par exemple, l'acetonitrile pur).

Prenez des mesures appropriées pour éviter les charges électrostatiques à proximité du système HPLC. Par exemple, s'assurez qu'il y a une humidité de l'air suffisante et une ventilation adéquate dans la laboratoire, portez des vêtements ou équipement de protection antistatique, évitez l'accumulation de bulles d'air dans les lignes de déchets et utilisez des réservoirs à déchets mis à la terre.

Utilisez uniquement des capillaires non-conducteurs pour diriger solvants au réservoir de déchets. Capillaires électriquement conducteur devrait être mis à la terre.

- **Inflammation spontanée des solvants**

N'utilisez aucun solvants avec une température d'auto-inflammabilité inférieure à 150° C. Si une fuite se produit, ces solvants peuvent s'auto-enflammer au contact d'une surface chaude.

- **Capillaires, connecteur capillaires, connexions ouvertes**

- ◆ Des capillaires, en particulier les capillaires non-métalliques, pourraient fendre ou glisser des connecteurs ou ne peuvent pas être vissés. Ceci peut en résulter aussi que des substances pourraient jaillir des connexions ouvertes.
- ◆ 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. Ceci peut causer des capillaires de fuite ou risquer des capillaires d'éclater. Ces acides peuvent cependant être utilisés dans le cadre de procédures de nettoyage, à condition que l'exposition soit brève.
- ◆ N'utilisez pas de capillaires écrasés, pliés, abimés ou endommagés.
- ◆ Les connecteurs capillaires pour pourrait être contaminé par des substances dangereuses ou des substances dangereuses pourrait sortir des connexions ouvertes.

- ◆ Certains capillaires des pompes RS, ainsi que des capillaires du système Viper, sont faits d'alliage de nickel-cobalt MP35N. Contact avec la peau peut provoquer une réaction chez les personnes qui sont sensibles au nickel/cobalt.
- ◆ Portez des lunettes de protection lorsque vous manipulez des capillaires en silice fondue (pendant l'installation, découpe, etc.).
- Pour raison de sécurité, ne soulevez pas l'instrument avec des réservoirs.
- Utilisez seulement des pièces de rechange originales et des accessoires autorisés par Thermo Fisher Scientific.
- Réglez toujours une limite de pression minimum pour le système HPLC. Ceci prévient les dommages résultant de fuites ou de long-terme fonctionnement à sec de la pompe. Activez la surveillance de niveau liquide pour des réservoirs (→ *manuel de pompe*).
- Lorsque vous soulevez ou l'instrument, tenez-le toujours par le dessous ou par les côtés de l'unité. Soulever l'instrument par la partie avant inférieure ou par le panneau avant peut endommager la porte.
- Ne placez aucun objet lourd sur la porte ouverte du panneau avant. Ceci pourrait endommager la porte.
- N'utilisez pas l'instrument de manière autre que celles décrites dans ce manuel.
- Conservez ce manuel à proximité de l'instrument pour pouvoir le consulter facilement.

1.3 Intended Use

For Research Use Only. Not for use in diagnostic procedures.

The device is designed to be operated only by qualified and authorized personnel. All users must know the hazards presented by the device and the used substances.

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 Installation and Maintenance

2.1 General Information

The solvent rack is a high-quality instrument designed for HPLC analysis as part of an UltiMate 3000 system. The rack is intended for use as a solvent organizer, for example, together with a LPG-3400 pump, providing a secure location for installation of the solvent reservoirs and, when mounted on top of the UltiMate 3000 system, saves valuable bench space.

All parts in the flow path are made of PEEK, FEP, stainless steel, or Titanium (biocompatible system only) to provide optimum resistance to the most commonly used HPLC solvents and buffer solutions.

2.2 Unpacking

All mechanical components of the solvent rack are carefully tested before the instrument is shipped from the factory. After unpacking, inspect the rack for any signs of mechanical damage that may have occurred during transit.

i **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 the packing material. They will provide excellent protection for the solvent rack in case of future transit. Shipping the rack in any other packaging automatically voids the product warranty. For more information, see the warranty statement in the terms of sale.

1. Place the shipping container on the floor. Remove the accessories kit and solvent bottles.
2. Grasp the solvent rack by the sides. Slowly and carefully, pull the rack out of the shipping container and place it on a stable surface.

⚠ Important: To prevent the rack from falling, always lift the unit by the sides. Do not lift the unit by the packaging material.

⚠ Important: Afin d'empêcher l'instrument de tomber, saisissez-la par les côtés. Ne soulevez l'instrument à l'aide du matériau d'emballage.

3. Remove the foam spacers, and then remove the polythene packaging.

2.3 Positioning the Solvent Rack in the UltiMate 3000 System

If the solvent rack is part of an UltiMate 3000 system, for example, for analytical HPLC applications, you should stack the individual modules, for example, as shown in Fig. 1. However, the arrangement of the system modules depends on the application.

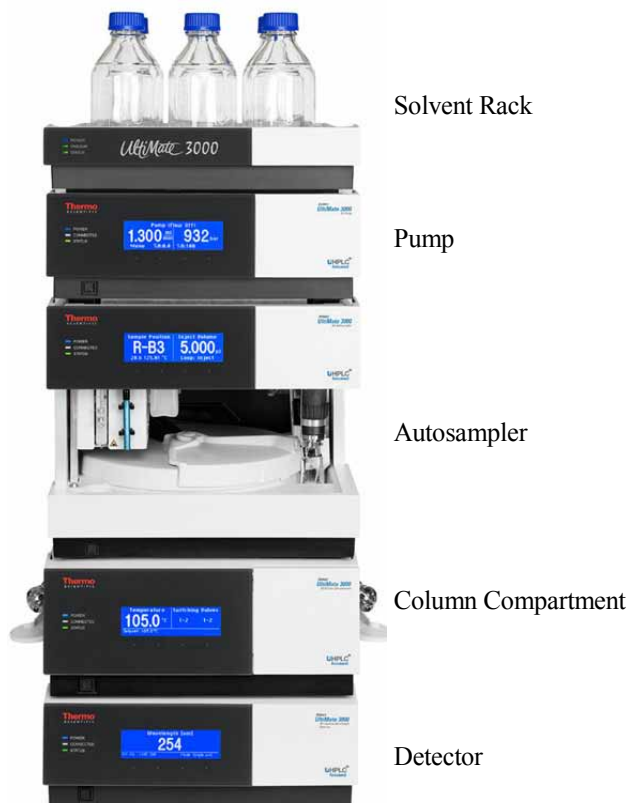


Fig. 1: Module arrangement for an UltiMate 3000 system (example)

If the solvent rack is part of an UltiMate 3000 RSLCnano or Proteomics MDLC system, the related system manual provides information about how to arrange the system modules and set up the systems for these applications.

2.4 Inside View

The front panel door tilts upward to provide access to the tubing guides in the solvent rack. The open door locks in its topmost position.

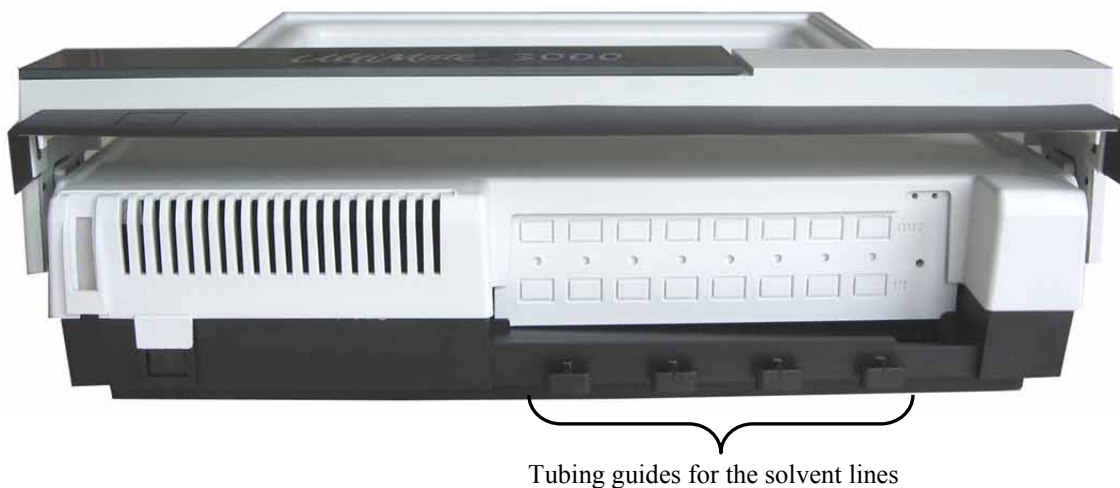


Fig. 2: SR-3000 with open front panel

- STOP Warning:** For safety reasons, you shall not lift the solvent rack when one or more solvent reservoirs are in the tray.
- STOP Avertissement:** Pour raison de sécurité, ne soulevez pas l'instrument avec des réservoirs.
- ! Important:** When lifting or moving the solvent rack, always lift by the bottom or sides of the unit. Lifting the solvent rack by the front panel may damage the front panel door.
- The open front panel door is not designed to carry weight. Therefore, you should not place any objects on the open door.
- ! Important:** Lorsque vous soulevez ou l'instrument, tenez-le toujours par le dessous ou par les côtés de l'unité. Soulever l'instrument par la partie avant inférieure ou par le panneau avant peut endommager la porte.
- Ne placez aucun objet sur la porte ouverte du panneau avant. Ceci peut endommager la porte.

2.5 Solvent Reservoirs

The solvent rack is shipped with solvent reservoirs and appropriate tubing, including filter holders and filter frits.

2.5.1 General Precautions

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

- Rinse the eluent reservoirs thoroughly before using them for the first time. Use high-purity solvent for rinsing.
- 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 *stainless steel* frits. If the UltiMate 3000 system includes a biocompatible pump, NCS-3500RS, or NCP-3200RS, replace the stainless frits with the frits from the accessories kit of the module. 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).

If the UltiMate 3000 system includes a semipreparative pump, use the filter holder and filter frit from the accessories kit of the pump.

- Locate the solvent reservoirs as close as possible to the pump. To avoid formation of gas bubbles in the reservoirs, 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. 1 (→ page 12). The tubing connecting the pump to the degasser should be as short as possible to avoid reformation of gas bubbles in the solvent.
- When replacing solvents, make sure that the solvents are miscible. Mix immiscible solvents with an intermediate solvent (for example, isopropanol) to replace them step-by-step.

2.5.2 Connecting the Solvent Reservoirs

1. Feed the solvent line through the retaining guide, which holds the tubing in place in the reservoir, and then into the open hole in the reservoir cap.
2. Verify that an appropriate frit is installed in the filter holder:

Pump	Filter Frit	Part No.
SD and SDN pumps	Stainless steel, porosity: 10 µm	6268.0110
BM pumps NCS-3500RS/NCP-3200RS	PEEK, porosity: 10 µm	6268.0117
RS, BX, and AB	Titanium, porosity: 10 µm	6268.0111

3. 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.
4. Place the complete assembly in the solvent reservoir.
5. Tighten the reservoir cap hand-tight. Press the retaining guide into the hole in the reservoir cap to ensure that the solvent line is hold in place inside the reservoir.



Fig. 3: Connecting the solvent lines to the reservoir

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

6. To connect the solvent lines to the pump, direct them through the opening in the rack.



Fig. 4: Openings in the solvent rack

7. Connect the solvent lines, for example, to the internal degasser of a LPG-3400 pump or directly to the pump inlet. Place the solvent lines in the tubing guides (→ Fig. 2, page 13), and then connect them to the pump.

2.6 Routine and Preventive Maintenance

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

Should you need to return an instrument, contact Thermo Fisher Scientific Service for Dionex HPLC Products. An RMA (Return Material Authorization) number is required in order to track your instrument. Always use the original packaging and observe the packing instructions when shipping the solvent rack. Shipping the rack in anything other than the original packaging will void 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.

3 Technical Information

Solvent reservoir capacity:	Eight 1-L reservoirs <i>or</i> four 2.5-L reservoirs <i>or</i> two 5-L reservoirs <i>or</i> five 1-L reservoirs and one 5-L reservoir <i>or</i> three 2.5L and two 1-L reservoirs <i>or</i> three 1-L and two 2.5-L reservoirs <i>or</i> sixteen 0.5L reservoirs
Wetted parts:	PEEK™, FEP, and ETFE/ECTFE Inline filter frit: SST, PEEK, or Titanium
Environmental conditions:	Range of use: Indoor use Temperature: 10 °C to 35°C Air humidity: max. 80% rel. humidity, non-condensing Pollution degree: 2
Dimensions (h x w x d):	10 x 42 x 51 cm (3.9 x 16.5 x 20 in)
Emission sound pressure level:	No emission
Weight:	3.0 kg (6.6 lbs) without reservoirs

Technical information: October 2012

All technical specifications are subject to change without notice.

4 Accessories, Consumables, and Spare Parts

4.1 Standard Accessories

Accessories for the solvent rack 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. The part numbers always refer to the packing unit. For more information contact the Thermo Fisher Scientific sales organization for Dionex HPLC Products.

The following accessories are shipped with the instrument (the list is subject to change without notice):

Description	Quantity in kit	Part No.
SR-3000 accessories, including:		
Solvent filter, including: Filter holder (top and bottom parts) and filter frit (stainless steel, 10µm)	4	Included in 6268.0115 Included in 6268.0110
Bottle cap for solvent reservoirs (including caps to close the holes in the bottle cap)	4	Included in 6270.0013
Cap (to close the holes in the bottle cap)	20	Included in 6000.0047
Retaining guide (solvent reservoir)	4	Included in 6000.0042
1-L solvent reservoir (with cap)	4	Included in 2270.0012

4.2 Consumables and Spare Parts

Description	Part No.
Filter holder (top and bottom parts) for solvent filter (pack of 6 filter holders) (to be used with all pumps except HPG-3200BX)	6268.0115
Filter holder (top and bottom parts) for solvent filter (pack of 2 filter holders) (to be used with all HPG-3200BX pumps)	6268.0116
Filter frit (stainless steel, 10µm) for solvent filter (pack of 10 frits) (to be used with SD and SDN pumps)	6268.0110
Filter frit (titanium, 10µm) for solvent filter (pack of 10 frits) (for example, to be used with RS, BX, and AB pumps)	6828.0111
Filter frit (PEEK, 10µm) for solvent filter (pack of 10 frits) (to be used with BM pumps and NCS-3500RS/NCP-3200RS)	6268.0117
Bottle cap for solvent reservoirs (including caps to close the holes in the bottle cap)	6270.0013
Cap to close the holes in the bottle cap (pack of 20 caps)	6000.0047
Retaining guide (solvent reservoir) (pack of 5 guides)	6000.0042
Solvent reservoir, 1-L reservoir with cap	2270.0012

