

Thermo

Surveyor MS Pump & Surveyor MS Pump Plus

Hardware Manual

60053-97105 Revision E December 2013



Thermo
SCIENTIFIC

© 2013 Thermo Fisher Scientific Inc. All rights reserved.

Xcalibur is a registered trademark and Surveyor is a trademark of Thermo Fisher Scientific Inc in the United States.

Super Flangeless is a trademark of Upchurch Scientific.

The following are registered trademarks in the United States and other countries: Windows is a registered trademark of Microsoft Corporation. Teflon is a registered trademark of E. I. du Pont de Nemours & Co.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history: Revision A, Apr 2005; Revision B, Apr 2006; Revision C, Sep 2006; Revision D, Jan 2009; Revision E, Dec 2013

This revision contains cosmetic changes only. The cosmetic changes include a new title page and chapter icon. In addition, the pagination might differ slightly from the last revision.

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

Regulatory Compliance

Thermo Fisher Scientific performs complete testing and evaluation of its products to ensure full compliance with applicable domestic and international regulations. When the system is delivered to you, it meets all pertinent electromagnetic compatibility (EMC) and safety standards as described in the next section or sections by product name.

Changes that you make to your system may void compliance with one or more of these EMC and safety standards. Changes to your system include replacing a part or adding components, options, or peripherals not specifically authorized and qualified by Thermo Fisher Scientific. To ensure continued compliance with EMC and safety standards, replacement parts and additional components, options, and peripherals must be ordered from Thermo Fisher Scientific or one of its authorized representatives.

Surveyor MS Pump Plus

EMC Directive 89/336/EEC amended by 92/31/EEC and 93/68/EEC

EMC compliance has been evaluated by Underwriters Laboratories Inc.

EN 55011	1998	EN 61000-4-3	2002
EN 61000-3-2	1995, A1; 1998, A2; 1998, A14; 2000	EN 61000-4-4	1995, A1; 2001, A2; 2001
IEC 61000-3-2	2000	EN 61000-4-5	1995, A1; 2001
EN 61000-3-3	1995	EN 61000-4-6	1996, A1; 2001
IEC 61000-3-3	1994	EN 61000-4-11	1994, A1; 2001
EN 61326-1	1997		
EN 61000-4-2	1995 A1; 1998 A2; 2001	CISPR 11	1999, A1; 1999, A2; 2002
FCC Class A, CFR 47 Part 15 Subpart B: 2003			

Low Voltage Safety Compliance

This device complies with Low Voltage Directive 73/23/EEC and harmonized standard EN 61010-1:2001.

FCC Compliance Statement

THIS DEVICE COMPLIES WITH PART 15 OF THE FCC RULES. OPERATION IS SUBJECT TO THE FOLLOWING TWO CONDITIONS: (1) THIS DEVICE MAY NOT CAUSE HARMFUL INTERFERENCE, AND (2) THIS DEVICE MUST ACCEPT ANY INTERFERENCE RECEIVED, INCLUDING INTERFERENCE THAT MAY CAUSE UNDESIRE OPERATION.



CAUTION Read and understand the various precautionary notes, signs, and symbols contained inside this manual pertaining to the safe use and operation of this product before using the device.

Notice on Lifting and Handling of Thermo Scientific Instruments

For your safety, and in compliance with international regulations, the physical handling of this Thermo Fisher Scientific instrument *requires a team effort* to lift and/or move the instrument. This instrument is too heavy and/or bulky for one person alone to handle safely.

Notice on the Proper Use of Thermo Scientific Instruments

In compliance with international regulations: This instrument must be used in the manner specified by Thermo Fisher Scientific to ensure protections provided by the instrument are not impaired. Deviations from specified instructions on the proper use of the instrument include changes to the system and part replacement. Accordingly, order replacement parts from Thermo Fisher Scientific or one of its authorized representatives.

Notice on the Susceptibility to Electromagnetic Transmissions

Your instrument is designed to work in a controlled electromagnetic environment. Do not use radio frequency transmitters, such as mobile phones, in close proximity to the instrument.

For manufacturing location, see the label on the instrument.

WEEE Compliance

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling or disposal companies in each European Union (EU) Member State, and these companies should dispose of or recycle this product. See www.thermoscientific.com/rohsweee for further information on Thermo Fisher Scientific's compliance with these Directives and the recyclers in your country.

WEEE Konformität

Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2002/96/EC erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet:








Thermo Fisher Scientific hat Vereinbarungen mit Verwertungs-/Entsorgungsfirmen in allen EU-Mitgliedsstaaten getroffen, damit dieses Produkt durch diese Firmen wiederverwertet oder entsorgt werden kann. Mehr Information über die Einhaltung dieser Anweisungen durch Thermo Fisher Scientific, über die Verwerter, und weitere Hinweise, die nützlich sind, um die Produkte zu identifizieren, die unter diese RoHS Anweisung fallen, finden sie unter www.thermoscientific.com/rohsweee.







Conformité DEEE

Ce produit doit être conforme à la directive européenne (2002/96/EC) des Déchets d'Équipements Électriques et Électroniques (DEEE). Il est marqué par le symbole suivant:



Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur www.thermoscientific.com/rohsweee.

CAUTION Symbol	CAUTION	VORSICHT	PRECAUCIÓN	MISE EN GARDE
	Risk electric shock: This instrument uses voltages that can cause electric shock and/or personal injury. Before servicing, shut down the instrument and disconnect it from line power. While operating the instrument, keep covers on. Do not remove the protective covers from the printed circuit board assemblies (PCBAs).	Stromschlaggefahr: Dieses Gerät arbeitet mit Spannungen, die Stromschläge und/oder Personenverletzungen verursachen können. Vor Wartungsarbeiten muss das Gerät abgeschaltet und vom Netz getrennt werden. Betreiben Sie das Gerät nicht mit abgenommenen Abdeckungen. Nehmen Sie die Schutzabdeckungen von Leiterplatten nicht ab.	Riesgo de descargas eléctricas: Este instrumento utiliza voltajes que pueden causar descargas eléctricas y/o lesiones personales. Antes de revisar o reparar el instrumento, apáguelo y desconéctelo de la red eléctrica. Mantenga colocadas las cubiertas mientras se utiliza el instrumento. No retire las cubiertas protectoras del circuito impreso completo (PCBA).	Risque de choc électrique : l'instrument utilise des tensions susceptibles de provoquer une électrocution et/ou des blessures corporelles. Il doit être arrêté et débranché de la source de courant avant toute intervention. Ne pas utiliser l'instrument sans ses couvercles. Ne pas enlever les capots de protection des cartes à circuit imprimé (PCBA).
	Chemical hazard: Wear gloves and other protective equipment, as appropriate, when handling toxic, carcinogenic, mutagenic, corrosive, or irritant chemicals. Use approved containers and proper procedures to dispose of waste oil and when handling wetted parts of the instrument.	Gefahr durch Chemikalien: Tragen Sie beim Umgang mit toxischen, karzinogenen, mutagenen, ätzenden oder reizenden Chemikalien Schutzhandschuhe und weitere geeignete Schutzausrüstung. Verwenden Sie bei der Entsorgung von verbrauchtem Öl und beim Umgang mit medienberührenden Komponenten die vorgeschriebenen Behälter, und wenden Sie ordnungsgemäße Verfahren an.	Peligro por sustancias químicas: Cuando manipule sustancias químicas, tóxicas, carcinogénicas, mutágenas, corrosivas o irritantes, utilice guantes y otro equipo de protección. Utilice siempre recipientes homologados y siga los procedimientos adecuados cuando deseche aceite residual o manipule partes mojadas del instrumento.	Danger lié aux produits chimiques : porter des gants et d'autres équipements de protection appropriés pour manipuler les produits chimiques toxiques, cancérogènes, mutagènes, corrosifs ou irritants. Utiliser des récipients homologués et des procédures adéquates pour la mise au rebut des huiles usagées et lors de la manipulation des pièces de l'instrument en contact avec l'eau.
	Hot surface: Before touching, allow any heated components to cool.	Heiße Oberflächen: Lassen Sie heiße Komponenten vor der Berührung abkühlen.	Superficies calientes: Antes de tocar los componentes calientes, espere a que se enfríen.	Surface chaude : laisser refroidir les composants chauffés avant toute manipulation.
	Flammable substances hazard: Use care when operating the system in the presence of flammable substances.	Gefahr durch entzündbare Substanzen: Beachten Sie die einschlägigen Vorsichtsmaßnahmen, wenn Sie das System in Gegenwart von entzündbaren Substanzen betreiben.	Peligro por sustancias inflamables: Tenga mucho cuidado cuando utilice el sistema cerca de sustancias inflamables.	Danger lié aux substances inflammables : agir avec précaution lors de l'utilisation du système en présence de substances inflammables.
	Risk of eye injury: Eye injury could occur from splattered chemicals, airborne particles, or sharp objects. (Sharp objects that customers might install in the instrument include fused-silica tubing, the autosampler needle, and so on.) Wear safety glasses when handling chemicals or servicing the instrument.	Augenverletzungsrisiko: Verspritzte Chemikalien, Schwebstoffpartikel oder scharfe Objekte können Augenverletzungen verursachen. (Scharfe Objekte, die Kunden möglicherweise im Gerät installieren, sind z. B. Quarzglas-Kapillaren, die Nadel des Autosamplers, usw.) Tragen Sie beim Umgang mit Chemikalien oder bei der Wartung des Gerätes eine Schutzbrille.	Riesgo de lesiones oculares: Las salpicaduras de sustancias químicas, las partículas flotantes en el aire y los objetos afilados pueden causar lesiones oculares. (Entre los objetos afilados que los clientes pueden instalar en el instrumento se encuentran tubos de sílice fundida, agujas del muestreador automático, etc.). Para manipular sustancias químicas o realizar tareas de mantenimiento, utilice gafas de seguridad.	Risque de lésion oculaire : les projections chimiques, les particules en suspension dans l'air et les objets tranchants peuvent entraîner des lésions oculaires. (Les objets tranchants pouvant être installés par les clients dans l'instrument comprennent les tubes en silice fondue, les aiguilles du passeur automatique, etc.). Porter des lunettes de protection lors de toute manipulation de produit chimique ou intervention sur l'instrument.
	General hazard: A hazard is present that is not included in the other categories. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual. When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.	Allgemeine Gefahr: Es besteht eine weitere Gefahr, die nicht in den vorstehenden Kategorien beschrieben ist. Dieses Symbol wird auch auf dem Gerät angebracht. Einzelheiten zu dieser Gefahr finden Sie in den Gerätehandbüchern. Wenn Sie sich über die Sicherheit eines Verfahrens im Unklaren sind, setzen Sie sich, bevor Sie fortfahren, mit dem technischen Support für Thermo Scientific San Jose Produkte in Verbindung.	Peligro general: Existen peligros que no se incluyen en las otras categorías. Este símbolo también aparece en el instrumento. Si desea obtener más información sobre estos peligros, consulte el manual del instrumento. En caso de duda sobre la seguridad de un procedimiento, póngase en contacto con el personal de servicio técnico de los productos Thermo Scientific San Jose.	Danger d'ordre général : indique la présence d'un risque n'appartenant pas aux catégories citées plus haut. Ce symbole figure également sur l'instrument. Pour plus de détails sur ce danger potentiel, se reporter au manuel de l'instrument. Si la sûreté d'une procédure est incertaine, contacter l'assistance technique pour les produits Thermo Scientific San Jose.

CAUTION Symbol	CAUTION	VORSICHT	PRECAUCIÓN	MISE EN GARDE
	Laser hazard: This instrument uses a laser that is capable of causing personal injury. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual.	Gefahr durch Laserstrahlen: Der in diesem Gerät verwendete Laser kann zu Verletzungen führen. Dieses Symbol wird auch auf dem Gerät angebracht. Einzelheiten zu dieser Gefahr finden Sie in den Gerätehandbüchern.	Peligro por láser: Este instrumento utiliza un láser que puede producir lesiones personales. Este símbolo también aparece en el instrumento. Si desea obtener más información sobre el peligro, consulte el manual del instrumento.	Danger lié au laser : l'instrument utilise un laser susceptible de provoquer des blessures corporelles. Ce symbole figure également sur l'instrument. Pour plus de détails sur ce danger potentiel, se reporter au manuel de l'instrument.
	Ultra violet light hazard: Do not look directly at the ultra-violet (UV) light or into the UV source. Exposure can cause eye damage. Wear UV eye protection.	Gefahr durch UV-Licht: Richten Sie Ihren Blick nicht direkt auf ultraviolettes Licht (UV-Licht) oder in die UV-Quelle. Dies kann zu Augenschäden führen. Tragen Sie eine UV-Schutzbrille.	Peligro por luz ultravioleta: No mire directamente a una luz ultravioleta (UV) ni a una fuente UV. La exposición puede causar daños oculares. Lleve protección ocular para UV.	Danger lié aux rayons ultraviolets : ne jamais regarder directement la lumière ultraviolette (UV) ou la source d'UV. Une exposition peut entraîner des lésions oculaires. Porter des protections oculaires anti-UV.
	Sharp object: Avoid physical contact with the object.	Scharfes Objekt: Vermeiden Sie den physischen Kontakt mit dem Objekt.	Objeto puntiagudo: Evite el contacto físico con el objeto.	Objet tranchant : éviter tout contact physique avec l'objet.
	Pinch point: Keep hands away from this area.	Quetschgefahr: Halten Sie Ihre Hände von diesem Bereich fern.	Puntos de pincamiento: Mantenga las manos apartadas de esta área.	Risque de pincement : éloigner les mains de cette zone.
	Heavy objects: Never lift or move the instrument by yourself; you can suffer personal injury or damage the equipment. For specific lifting instructions, refer to the instrument manual.	Schweres Objekt: Bewegen und heben Sie das Gerät niemals allein an; dies kann zu Verletzungen oder zur Beschädigung des Geräts führen. Spezifische Anweisungen zum Anheben finden Sie im Gerätehandbuch.	Objeto pesado: Nunca levante ni mueva el instrumento por su cuenta, podría sufrir lesiones personales o dañar el equipo. Para obtener instrucciones específicas sobre levantamiento, consulte el manual del instrumento.	Objet lourd : ne jamais soulever ou déplacer l'instrument seul sous peine de blessure corporelle ou d'endommagement de l'instrument. Pour obtenir des instructions de levage spécifiques, se reporter au manuel de l'instrument.
	Trip obstacle: Be aware of cords, hoses, or other objects located on the floor.	Stolpergefahr: Achten Sie auf Kabel, Schläuche und andere Objekte auf dem Fußboden.	Tropiezo con obstáculos: Tenga en cuenta los cables, mangueras u otros objetos colocados en el suelo.	Risque de trébuchement : faire attention aux câbles, tuyaux et autres objets situés sur le sol.
	When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.	Wenn Sie sich über die Sicherheit eines Verfahrens im unklaren sind, setzen Sie sich, bevor Sie fortfahren, mit Ihrer lokalen technischen Unterstützungsorganisation für Thermo Scientific San Jose Produkte in Verbindung.	En caso de duda sobre la seguridad de un procedimiento, póngase en contacto con el personal de servicio técnico de los productos Thermo Scientific San Jose.	Si la sûreté d'une procédure est incertaine, contacter l'assistance technique pour les produits Thermo Scientific San Jose.

CAUTION Symbol	CAUTION	警告	危険警告
	<p>Risk electric shock: This instrument uses voltages that can cause electric shock and/or personal injury. Before servicing, shut down the instrument and disconnect it from line power. While operating the instrument, keep covers on. Do not remove the protective covers from the printed circuit board assemblies (PCBAs).</p>	<p>感電の危険性: この機器では、感電および/または身体傷害を引き起こすおそれのある電圧を使用しています。整備点検の前には、機器の電源を切り、電源コードを抜いてください。機器の作動中は、カバーを付けたままにしてください。プリント基板アセンブリ (PCBA) から保護カバーを取り外さないでください。</p>	<p>触电危险: 本仪器所用电压可能导致电击或人身伤害。进行维修服务前, 务必关闭仪器电源并断开其电源连接。操作此仪器时, 不要卸下顶盖。勿卸下印刷电路板组件 (PCBA) 的保护盖。</p>
	<p>Chemical hazard: Wear gloves and other protective equipment, as appropriate, when handling toxic, carcinogenic, mutagenic, corrosive, or irritant chemicals. Use approved containers and proper procedures to dispose of waste oil and when handling wetted parts of the instrument.</p>	<p>化学的危険性: 毒性、発癌性、変異原性、腐食性、または刺激性のある化学薬品を取り扱うときは、必要に応じて手袋などの保護具を着用します。廃油を処分したり、機器の接液部品を取り扱うときは、認可された容器を使用し、適切な手順に従います。</p>	<p>化学品危险: 当处理毒性、致癌性、致突变性、腐蚀性或者刺激性化学品时, 佩戴手套和其他保护性设备。当处理浸湿的仪器部件以及废油时, 使用认可的容器和合适的步骤。</p>
	<p>Hot surface: Before touching, allow any heated components to cool.</p>	<p>高温面: 触れる前に、加熱した部品を冷ましてください。</p>	<p>热表面: 待高温部件冷却之后再行维修。</p>
	<p>Flammable substances hazard: Use care when operating the system in the presence of flammable substances.</p>	<p>可燃性物質の危険性: 可燃性物質があるところでシステムを作動させる場合は十分注意してください。</p>	<p>易燃物危险: 在有易燃物质的场地操作该系统时, 务必小心谨慎。</p>
	<p>Risk of eye injury: Eye injury could occur from splattered chemicals, airborne particles, or sharp objects. (Sharp objects that customers might install in the instrument include fused-silica tubing, the autosampler needle, and so on.) Wear safety glasses when handling chemicals or servicing the instrument.</p>	<p>眼外傷の危険性: 飛散した化学薬品、浮遊粒子、または鋭利な物体によって眼外傷を負うおそれがあります (機器に取り付けられる可能性がある鋭利な物体は、ヒューズドシリカ、オートサンプラーニードルなどです)。化学薬品を取り扱ったり、機器を整備点検するときは、保護メガネを着用します。</p>	<p>眼睛伤害风险: 眼睛受伤可能源自飞溅的化学品、空气中的颗粒, 或者锋利的物体。(安装在仪器内的锋利物体包括熔融石英管、自动进样器的进样针等。) 处理化学品或对仪器进行维修服务时, 务必戴上防护眼镜。</p>
	<p>General hazard: A hazard is present that is not included in the other categories. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual. When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.</p>	<p>一般的な危険性: それぞれのカテゴリーに当てはまらない危険があります。この標識記号は機器にも表示されています。この危険の詳細については、機器のマニュアルを参照してください。手順の安全性にご不明な点がある場合は、Thermo Scientific San Jose 製品のテクニカルサポートまでお問い合わせください。</p>	<p>普通危険: 未归入其他类别的危险。此符号也会在仪器上出现。有关此危险的详细信息, 参阅适当的仪器手册。若对任何步骤的安全事项有疑问, 联系 Thermo Scientific San Jose 产品的技术支持中心。</p>

CAUTION Symbol	CAUTION	警告	危険警告
	Laser hazard: This instrument uses a laser that is capable of causing personal injury. This symbol also appears on the instrument. For details about the hazard, refer to the instrument manual.	レーザー光線の危険性: この機器では、身体傷害を引き起こすおそれのあるレーザーを使用しています。この標識記号は機器にも表示されています。この危険の詳細については、機器のマニュアルを参照してください。	激光危险: 本仪器所用激光会导致人身伤害。此符号也会在仪器上出现。有关此危险的详细信息, 参阅适当的仪器手册。
	Ultra violet light hazard: Do not look directly at the ultra-violet (UV) light or into the UV source. Exposure can cause eye damage. Wear UV eye protection.	紫外光の危険性: 紫外(UV)光またはUV光源を直接見ないでください。照射によって眼損傷を引き起こすおそれがあります。UV保護メガネを着用します。	紫外光危险: 不要直视紫外(UV)光或者紫外光源。直视可能导致眼睛伤害。佩戴紫外线防护眼镜。
	Sharp object: Avoid physical contact with the object.	鋭利な物体: 物体との身体的接触を避けてください。	锋利物体: 避免直接接触锋利的物体。
	Pinch point: Keep hands away from this area.	ピンチポイント: この部分には手を挟まれないようにしてください。	夹点: 勿将手放在此部位。
	Heavy objects: Never lift or move the instrument by yourself; you can suffer personal injury or damage the equipment. For specific lifting instructions, refer to the instrument manual.	重量物: 1人で機器を持ち上げたり移動しないでください。身体傷害を負ったり、機器を損傷するおそれがあります。具体的な持ち上げ方法については、機器のマニュアルを参照してください。	重物: 切勿独自提起或移动本仪器; 可能遭受人身伤害或损坏仪器。有关具体的提起说明, 参阅仪器手册。
	Trip obstacle: Be aware of cords, hoses, or other objects located on the floor.	作業の障害物: 床にあるコード、ホース、その他の物体に注意してください。	绊倒危险: 注意地面上的线、管或其他物品。
	When the safety of a procedure is questionable, contact Technical Support for Thermo Scientific San Jose products.	手順の安全性にご不明な点がある場合は、Thermo Scientific San Jose 製品のテクニカルサポートまでお問い合わせください。	如对安全程序有疑问, 联系 Thermo Scientific San Jose 产品的技术支持中心。

Contents

	Preface	xiii
	Related Documentation	xiii
	Cautions and Special Notices	xiv
	Good Laboratory Practices	xiv
	Contacting Us	xvi
Chapter 1	Introduction	1
	Functional Description	2
	Vacuum Degasser Assembly	3
	Liquid Displacement Assembly	3
	Pulse Dampening Assembly	4
	Status LEDs	6
	Specifications	7
Chapter 2	Installation	9
	Installation Checklist	10
	Unpacking and Inspecting the Instrument	10
	Making the Initial Instrument Preparations	11
	Making the Initial Back Panel Connections	11
	Serial Communication Cable	12
	System Interconnect Cable	12
	Connecting the Low-Pressure Solvent Lines	15
	Step 1: Assembling the Solvent Reservoir Bottles	16
	Step 2: Installing the Super Flangeless Fittings	17
	Connecting the Pump to the Surveyor Autosampler Plus	19
	Powering On for the First Time	21
	Configuring the Data System PC COM Port	21
	Preparing the Pump for Operation	22
	Specifying the Instrument Configuration Options	22
	Purging Air from the Solvent Lines	25
	Priming the Pulse Dampener	26

Chapter 3	Diagnostics	29
	Monitoring Pump Pulsation	29
	Testing the Pump Proportioning	31
	Determining the Delay Volume of the Pump	33
Chapter 4	Maintenance	35
	Flushing the Piston Guide Bushings	36
	Purging Air from the Piston Chambers	37
	Cleaning the LDA Components without Disassembling the LDA	38
	Servicing the LDA	39
	Removing the LDA from the Pump	40
	Disassembling the LDA	43
	Cleaning the LDA Components	47
	Replacing the Primary Piston Seals	49
	Reassembling the LDA	51
	Reinstalling the LDA	59
	Servicing the Proportioning Valves	62
	Replacing the Solvent Reservoir Filter	63
Chapter 5	Troubleshooting	65
	Troubleshooting Tips	65
	General System Troubleshooting	66
	Surveyor MS Pump Troubleshooting Guide	68
	Pump Pulsation	70
	Normal Pressure Pulsation	70
	Abnormal Pressure Pulsation	71
	Baseline Noise at Low Pressures (under 40 Bar)	71
	Insufficient Solvent Flow from the Pump Outlet	72
	Pump Cavitation	72
	Blocked Proportioning Valve	73
	Blocked Check Valve	73
	Irreproducible Retention Times	73
Chapter 6	Accessories and Replacement Parts	75
	Liquid Displacement Assembly (LDA) Parts	76
	MS Pump Accessory Kit	78
	Solvent Interconnect Kit	78
	Other Parts and Assemblies	79
	Index	81

Preface

The Surveyor MS Pump and Surveyor MS Pump Plus are members of the Surveyor Plus™ family of LC instruments.

This *Surveyor MS Pump and Surveyor MS Pump Plus Hardware Manual* provides you with information on how to setup, maintain, and troubleshoot the Surveyor MS Pump and the Surveyor MS Pump Plus. A list of accessories and spare parts is also included. The Surveyor MS Pump and the Surveyor MS Pump Plus are operated from a computer running the Xcalibur™ data system.

Contents

- [Related Documentation](#)
- [Cautions and Special Notices](#)
- [Contacting Us](#)

Related Documentation

In addition to this guide, Thermo Fisher Scientific provides the following documents for the Surveyor MS Pump and the Surveyor MS Pump Plus:

- *Surveyor Plus Preinstallation Requirements Guide*
- *Surveyor Plus Getting Connected Guide*
- *Surveyor Plus User Guide for LC Devices*

Cautions and Special Notices

Make sure you follow the cautions and special notices presented in this guide. Cautions and special notices appear in boxes; those concerning safety or possible system damage also have corresponding caution symbols.

This guide uses the following types of cautions and special notices.



CAUTION Highlights hazards to humans, property, or the environment. Each CAUTION notice is accompanied by an appropriate CAUTION symbol.

IMPORTANT Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the system.

Note Highlights information of general interest.

Tip Highlights helpful information that can make a task easier.

Good Laboratory Practices

To obtain optimal performance from your LC system and to prevent personal injury or injury to the environment, do the following:

- Keep good records.
- Read the manufacturers' Material Safety Data Sheets (MSDSs) for the chemicals being used in your laboratory.
- Remove particulate matter from your samples before you inject them into the liquid chromatograph.
- Use HPLC grade solvents.
- Connect the drainage tubes from the pump, autosampler, and detector to an appropriate waste receptacle. Dispose of solvents as specified by local regulations.

Keep Good Records

To help identify and isolate problems with either your equipment or your methodology, we recommend that you keep good records of all system conditions (for example, % RSDs on retention times and peak areas, peak shape and resolution). At a minimum, keep a chromatogram of a typical sample and standard mixture, well documented with system

conditions, for future reference. Careful comparison of retention times, peak shapes, peak sensitivity, and baseline noise can provide valuable clues to identifying and solving future problems.

Chemical Toxicity

Although the large volume of toxic and flammable solvents used and stored in laboratories can be quite dangerous, do not ignore the potential hazards posed by your samples. Take special care to read and follow all precautions that ensure proper ventilation, storage, handling, and disposal of both solvents and samples. Become familiar with the toxicity data and potential hazards associated with all chemicals by referring to the manufacturers' Material Safety Data Sheets.

Sample Preparation

Always consider the solubility of your sample in the solvent/mobile phase. Sample precipitation can plug the column, tubing or flow cell causing flow restriction. This obstruction can result in irreparable damage to the system. Particulate matter can be avoided by filtering the samples through 0.45 or 0.2 micron (or less) filters.

Solvent Requirements

Many chemical manufacturers provide a line of high-purity or HPLC-grade reagents that are free of chemical impurities. Routine filtration of all solvents or eluents through a 0.45 or 0.2 micron (or less) fluorocarbon filter before placing them in the solvent reservoir significantly prolongs the life and effectiveness of the inlet filters, check valves and seals, injector, and column. Typically, HPLC-grade solvents do not require filtration.

Choose a mobile phase that is compatible with the sample and column you have selected for your separation. Remember that some solvents are corrosive to stainless steel.

Solvent Disposal

Make sure you have a solvent waste container or other kind of drain system available at or below the benchtop level. Most solvents have special disposal requirements and should not be disposed of directly down a drain. Follow all governmental regulations when disposing of any chemical.

High-pressure Systems and Leaks

LC systems operate at high pressures. Because liquids are not highly compressible they do not store much energy. Accordingly, there is little immediate danger from the high pressures in an LC system. However, if a leak occurs, it should be corrected as soon as possible. Always wear eye and skin protection when operating or maintaining an LC system. Always shut down the system and return it to atmospheric pressure before attempting any maintenance.

Contacting Us

There are several ways to contact Thermo Fisher Scientific for the information you need.

❖ **To contact Technical Support**

Phone	800-532-4752
Fax	561-688-8736
E-mail	us.techsupport.analyze@thermofisher.com
Knowledge base	www.thermokb.com

Find software updates and utilities to download at mssupport.thermo.com.

❖ **To contact Customer Service for ordering information**

Phone	800-532-4752
Fax	561-688-8731
E-mail	us.customer-support.analyze@thermofisher.com

❖ **To get local contact information for sales or service**

Go to <http://www.thermoscientific.com/en/support-landing/support.html>.

❖ **To copy manuals from the Internet**

Go to mssupport.thermo.com, agree to the Terms and Conditions, and then click **Customer Manuals** in the left margin of the window.

❖ **To suggest changes to the documentation or to the Help**

- Fill out a reader survey online at www.surveymonkey.com/s/PQM6P62.
- Send an e-mail message to the Technical Publications Editor at techpubs-lcms@thermofisher.com.

Introduction

The Surveyor MS Pump and Surveyor MS Pump Plus are members of the Surveyor Plus™ family of LC instruments.

This chapter provides an introduction to the Surveyor MS Pump and Surveyor MS Pump Plus (see [Figure 1](#)).

Thermo Fisher Scientific no longer ships the Surveyor MS Pump, but this legacy version of the pump can be controlled with the new instrument control software provided on the LC Devices software CD. The information provided in this manual applies to both the Surveyor MS Pump and the Surveyor MS Pump Plus.

Contents

- [Functional Description](#)
- [Specifications](#)

Figure 1. Surveyor MS Pump Plus



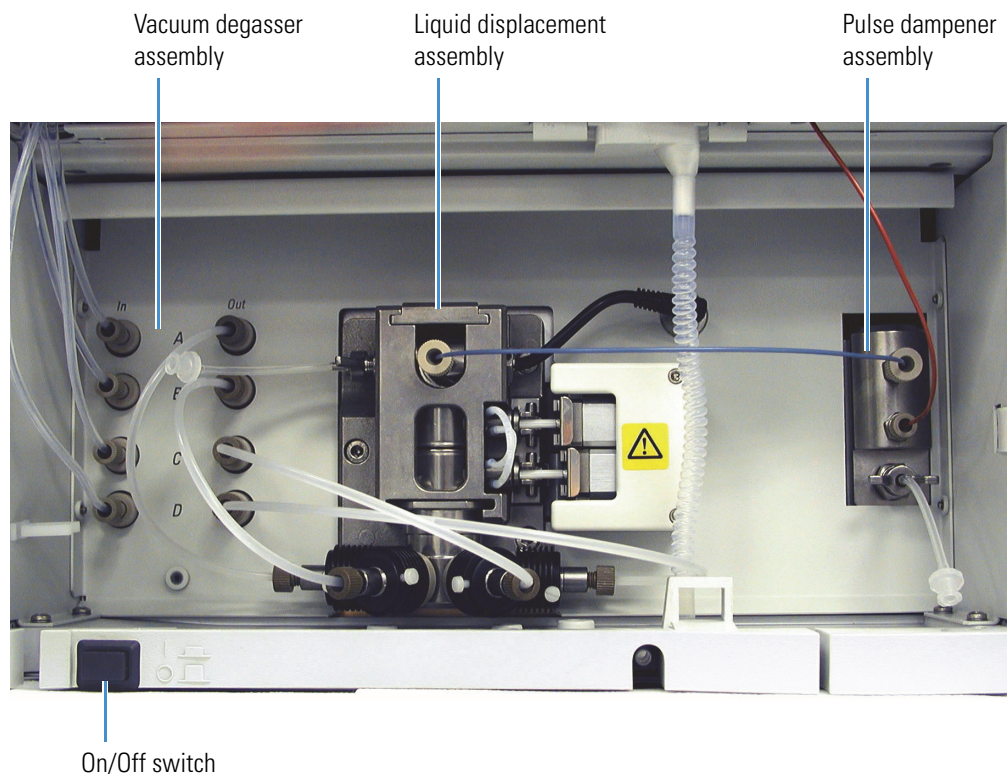
Functional Description

The Surveyor MS Pump Plus is a quaternary, low-pressure mixing pump with built-in solvent degassing and pulse dampening systems. The pumping system provides optimum performance in the lower flow rate ranges needed for mass spectrometry. You can run precise gradients from 25 to 2000 $\mu\text{L}/\text{min}$. An extremely low gradient delay volume of 80 μL ensures minimum system cycle times. The integral vacuum degasser offers superb solvent degassing efficiency with less than 500 μL of volume, and the pulse dampener produces stable flow rates with only 3 μL of delay volume.

The Surveyor MS Pump Plus is a bench-top unit for inclusion in the Surveyor high performance liquid chromatograph (HPLC). It is remotely controlled by way of an RS232 serial communication link from a computer using Microsoft Windows NT-based Xcalibur software (version 1.3 or higher). The only manual control on the pump is a power switch.

The Surveyor MS Pump Plus (see [Figure 2](#)) has three major components: the vacuum degasser assembly, the liquid displacement assembly (LDA), and the pulse dampening assembly. In addition, the MS pump contains status LEDs, a low voltage power supply, and several printed circuit boards (PCBs).

Figure 2. Surveyor MS Pump Plus with the doors open, showing the major components



Vacuum Degasser Assembly

The Surveyor MS Pump Plus contains a built-in vacuum membrane degasser for the purpose of removing dissolved gases from the mobile phase. Dissolved gases can cause many problems in HPLC and must be kept to an absolute minimum for best performance. The MS pump solvent degassing system consists of four independent channels that are maintained at a constant vacuum of approximately 686 mm (27 in.) Hg. Each channel has a volume of less than 500 μL , which provides superior degassing efficiency capable of limiting dissolved oxygen to a level of only 0.8 ppm. The system's vacuum pump operates continuously, which eliminates the excess baseline noise and drift caused by vacuum pump cycling.

Liquid Displacement Assembly

The frame of the liquid displacement assembly (LDA) consists of two components: the LDA body and the inlet module. The LDA body contains the inlet and outlet check valves, two TZP-ceramic pistons and their seals, an upper and lower piston chamber, a pressure transducer, and various PTFE seal rings (see [Figure 75 on page 77](#)). The inlet module contains a built-in low volume mixer. Four high-precision proportioning valves are connected to the inlet module. And four actuating coils are connected to the ends of the proportioning valves. All of these components are integrated into a single mechanical-hydraulic assembly, which minimizes the pump's gradient delay volume to only 80 μL .

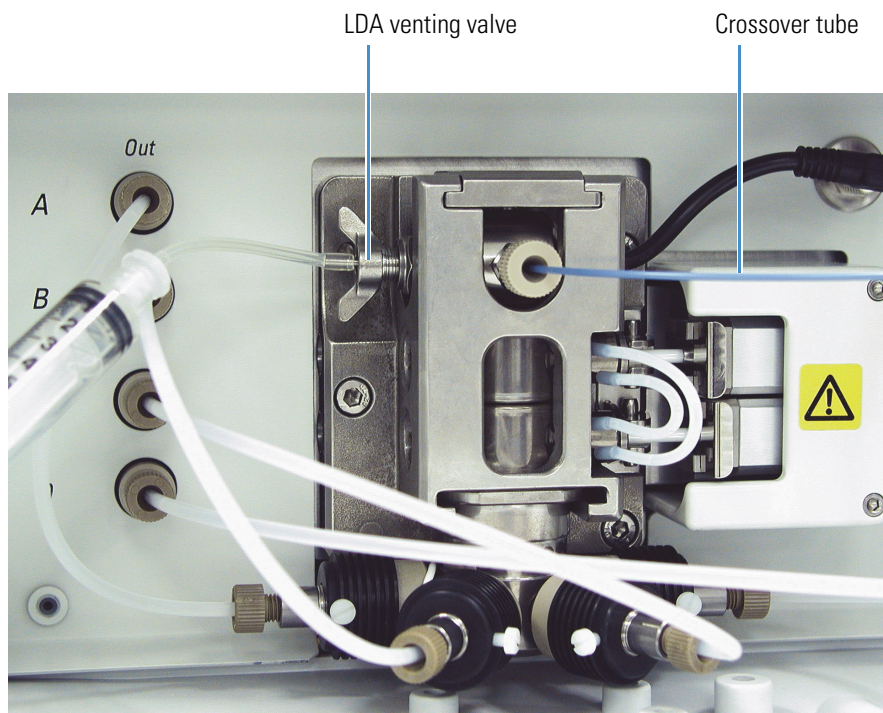
As the lower piston withdraws, mobile phase from the proportioning valve assembly passes through the inlet check valve and fills the lower piston chamber (see [Figure 75 on page 77](#)). As the lower piston returns, the inlet check valve closes and the mobile phase is pushed through the outlet check valve and into the upper piston chamber. Both check valves are low volume and incorporate a dual ball and seat design, which ensures a positive seal with minimal delay volume.

Next, the upper piston moves into the upper piston chamber, closing the outlet check valve and forcing the mobile phase into the pressure transducer housing. On the next piston stroke the mobile phase is pushed out of the LDA and into the crossover tube that is attached to the pulse dampening assembly.

The proportioning valves of the Surveyor MS Pump Plus incorporate a finely machined ball and seat design similar to a check valve. This ball and seat design controls the flow of solvents to a higher degree of precision and accuracy than the diaphragm design used by most gradient pumps. In addition, the ruby ball and sapphire seat are virtually wear free as opposed to the polymeric membrane of a diaphragm valve, which becomes more brittle and less responsive as it ages.

There is a venting valve on the left side of the LDA. When the venting valve is open, the mobile phase exits out the left end of the pressure sensor assembly. When the venting valve is closed, the mobile phase is directed out of the front of the pressure sensor assembly and into the crossover tube. [Figure 3](#) shows a pump with the venting valve in the open position for purging.

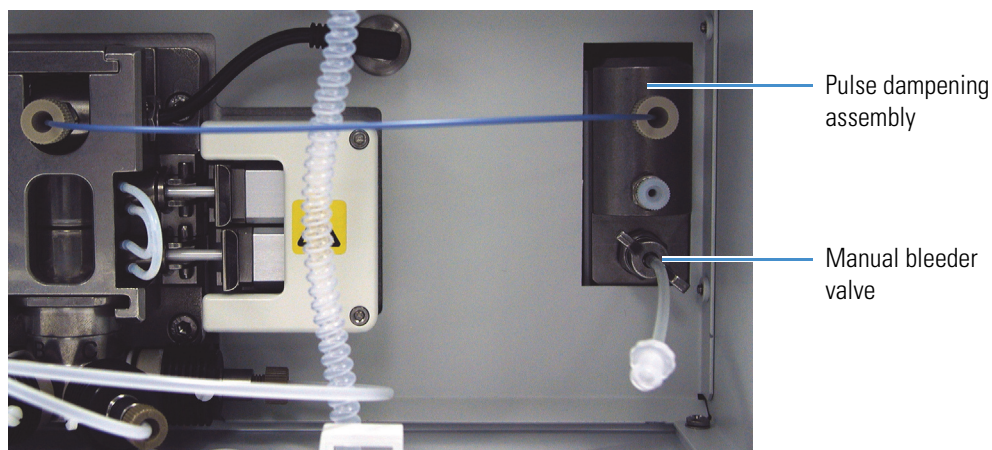
Figure 3. Purging air from the Surveyor MS Pump Plus



Pulse Dampening Assembly

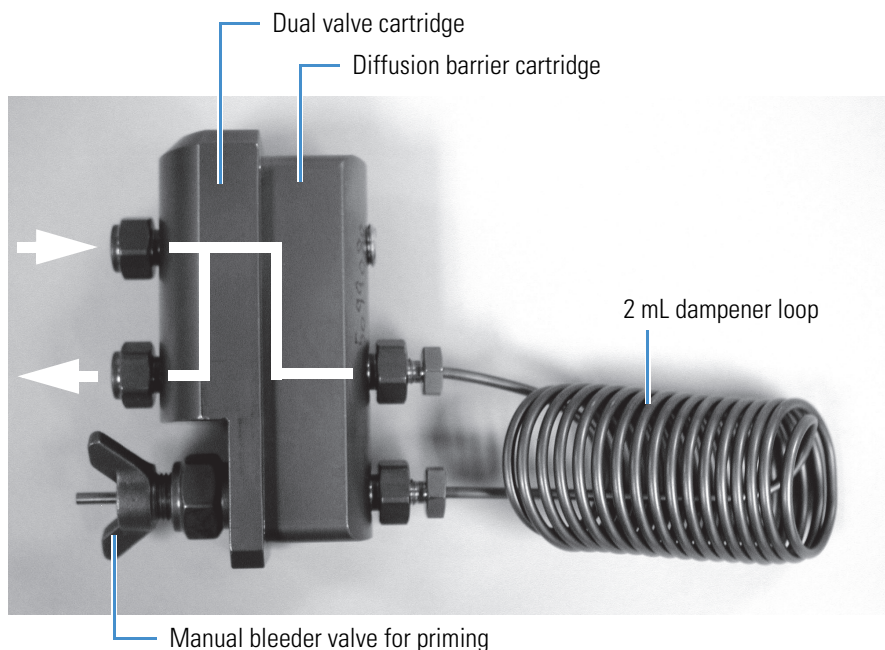
The pulse dampening assembly is located on the right side of the Surveyor MS Pump Plus. See [Figure 4](#).

Figure 4. Front view of the pulse dampening assembly



The pulse dampening assembly, built into the Surveyor MS Pump Plus, consists of a low volume T-connector through which the mobile phase passes. Attached to the side leg of the T-connector is a permeable isolation membrane made of sintered Teflon followed by a 2 mL loop of stainless steel tubing. See Figure 5.

Figure 5. Pulse dampening assembly



The loop is terminated with a priming valve. When the priming valve is open, the dampening loop can be flushed or filled with an appropriate solvent such as methanol or isopropanol. When the priming valve is closed, the diffusion barrier cartridge is shutoff from the flow path and the dampening loop absorbs pump pulsations. Even at elevated pressures, the pump shows only minimal pulsation when the pulse dampener is used. And because the 2 mL loop is shut off from the flow path, the low volume T-connector of the pulse dampener adds only 3 μL of delay volume to the system.

Because the membrane between the T-connector and the 2 mL loop is permeable, pumping buffered mobile phases leads to a slow buildup of electrolytes in the loop. To prevent corrosion in this situation, flush and prime the pulse dampener periodically.



CAUTION Do not open the manual bleeder valve during normal operation of the pump. If you open the valve while you are pumping an acidic or buffered mobile phase, prime the pulse dampener as soon as you notice the error. Because the pulse dampener is composed of two dissimilar metals, titanium and stainless steel, leaving acids or buffered solutions inside it causes galvanic corrosion and increased baseline noise.



CAUTION Do not confuse the manual bleeder valve with the venting valve on the left side of the LDA. The manual bleeder valve **is not** a purge valve for quickly drawing solvents through the solvent lines.

Status LEDs

Four status LEDs – labeled Power, Comm, Run, and Degas – are located on the right door of the pump. See [Figure 6](#). The states of these LEDs are described in [Table 1](#).

Figure 6. Front panel LEDs



Table 1. Status states for the Surveyor MS Pump Plus

LED	State	Meaning
Power	Amber	The MS Pump is switched On and is receiving power, but has not yet passed its Power On Self Test.
	Green	The MS Pump is switched On and is receiving power, and has passed its Power On Self Test.
Comm	Amber	The MS pump is not communicating with the data system PC
	Green	Communication to the data system PC has been established
	Flashing green	A pump program is downloading from the data system PC.
Run	Green	The MS pump is ready for a run.
	Flashing green	The MS pump is in a run.
Degas	Amber	The degas unit is building vacuum.
	Flashing amber	A failure, such as a loss of vacuum, has occurred.
	Green	Sufficient vacuum has developed to perform chromatography.

Specifications

The specifications for the Surveyor MS Pump Plus are listed below.

Minimum Flow Rate:	0.2 $\mu\text{L}/\text{min}$ isocratic, 25 $\mu\text{L}/\text{min}$ gradient
Maximum Flow Rate:	2000 $\mu\text{L}/\text{min}$
Flow Rate Resolution:	0.1 $\mu\text{L}/\text{min}$
Pressure Range:	0 to 400 bar (0 to 5800 psi)
Operating Range:	40 to 400 bar (588 to 5800 psi)
Pressure Resolution:	0.01 bar (0.15 psi)
Solvent Capacity:	Quaternary
Composition Accuracy:	$\pm 1\%$ at a flow rate of 200 $\mu\text{L}/\text{min}$
Piston Volume:	24 μL
Delay Volume:	80 μL
Degasser Type:	Vacuum membrane
Degasser Capacity:	4 channels
Degasser Volume:	< 500 μL per channel
Pulse Dampener:	3 μL , dynamic
Wetted Parts:	316 stainless steel, titanium, PEEK, sapphire, TZP-ceramic, FEP, GFP, ruby
Remote Controls:	Start, RS232 serial interface for Xcalibur software
Dimensions:	18 cm \times 36 cm \times 47 cm ($h \times w \times d$)
Weight:	12 kg
Power Requirements:	100/120 V ac or 220/240 V ac, 50 to 60 Hz, 500 VA max
Operating Temp:	+ 10 $^{\circ}\text{C}$ to + 40 $^{\circ}\text{C}$
Storage Temp:	- 40 $^{\circ}\text{C}$ to + 70 $^{\circ}\text{C}$
Operating Humidity:	5 to 95% relative, non-condensing

Installation

This chapter provides instructions for the initial installation of your Surveyor MS Pump Plus, including connection to other modules in the system. The Installation Checklist on the back of this page is an abbreviated version of the chapter and can be used as a quick reference of how to conduct a successful installation. Make a copy of the checklist and fill it out when the installation is complete. Include the completed checklist in your maintenance records.

Note Perform the installation in the sequence presented on the Installation Checklist and detailed in this chapter.

Contents

- [Installation Checklist](#)
- [Unpacking and Inspecting the Instrument](#)
- [Making the Initial Instrument Preparations](#)
- [Making the Initial Back Panel Connections](#)
- [Connecting the Low-Pressure Solvent Lines](#)
- [Connecting the Pump to the Surveyor Autosampler Plus](#)
- [Powering On for the First Time](#)
- [Preparing the Pump for Operation](#)

Installation Checklist

This topic provides a brief summary of the steps that must be completed for the proper installation of your Surveyor MS Pump Plus:

- Unpack and inspect your instrument (“[Unpacking and Inspecting the Instrument](#)” on [page 10](#)).
- Read the safety precautions.
- Position the pump appropriately (“[Making the Initial Instrument Preparations](#)” on [page 11](#)).
- Make the back panel connections (“[Making the Initial Back Panel Connections](#)” on [page 11](#)).
- Connect the solvent lines (“[Connecting the Low-Pressure Solvent Lines](#)” on [page 15](#)).
- Connect the power cord and turn on the instrument (“[Powering On for the First Time](#)” on [page 21](#)).
- Prime the pulse dampener (“[Priming the Pulse Dampener](#)” on [page 26](#)).

This Surveyor MS Pump Plus was installed by:

(Name)

(Date)

Unpacking and Inspecting the Instrument

Carefully remove the pump from the shipping container and inspect both the pump and packing for any signs of damage. If you find any damage, save the shipping materials and immediately contact the shipping company.

The shipping container should contain the pump, an accessory kit, a power cable, a serial communications cable with DB-9 pin connectors, and this manual. The accessory kit contains two 1/16 in. PEEK fittings, a 10 mL Luer-tipped syringe, a 4 mm balldriver with a screwdriver style handle, and a 4 mm Allen wrench.

Carefully check to make sure you received all the items listed on the packing list. If any items are missing, contact your Thermo Fisher Scientific service representative immediately.

Making the Initial Instrument Preparations

Place the pump on a laboratory bench as close as possible to a designated electrical outlet. Be sure to place the system in a **draft-free** location away from an open window, air conditioner vents, or other circulating air source. A stable room temperature is necessary for applications requiring maximum detection sensitivity. The area should also be free from dust, moisture, direct sunlight, strong electromagnetic fields, and physical vibrations. Allow at least 15 cm (6 in.) of clear space between the back panel of the pump and any wall or obstruction. This provides access to the back-panel connectors and a free flow of cooling air.



CAUTION Never connect the pump to, or operate the pump with, an electrical line source that has power drops or fluctuations of 10% below the nominal rated line voltage.

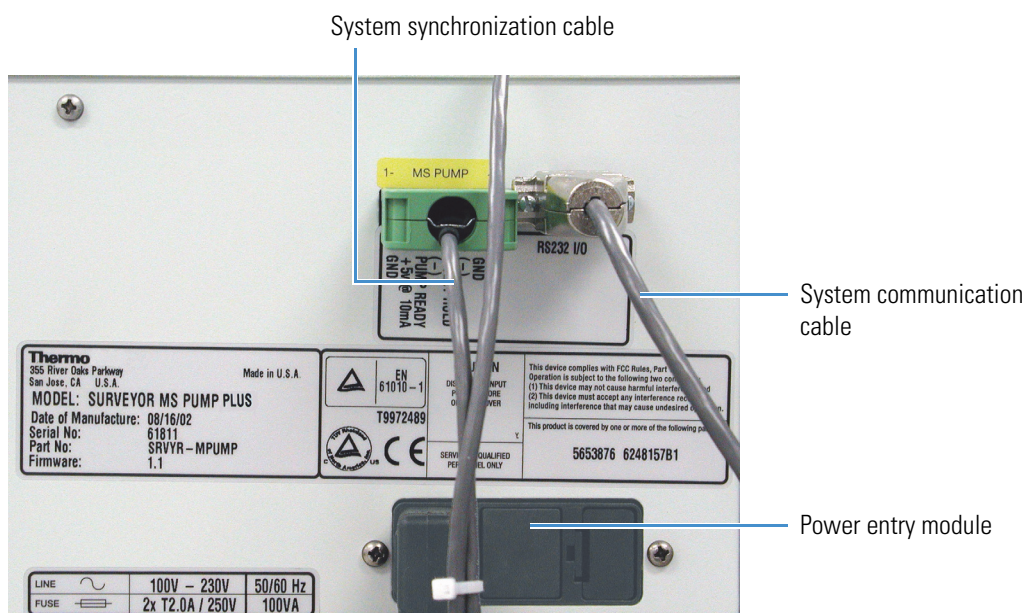


CAUTION Do not connect the power cord or turn on the pump yet. Wait until the [Powering On for the First Time](#) section of this chapter.

Making the Initial Back Panel Connections

In addition to connecting the MS pump to line power, operation of the Surveyor MS Pump Plus requires connecting two cables to the back panel of the pump. See [Figure 7](#). One is a serial communication cable that is used to download analytical methods to, and upload pressure information from, the pump. The other is a system synchronization cable that is used to coordinate the operation of the pump with the other Surveyor modules.

Figure 7. Back panel connections



2 Installation

Making the Initial Back Panel Connections

Serial Communication Cable

The serial communication cable is a standard RS232 cable, approximately 2 meters long and terminated at each end with a DB-9 pin connector. The serial communication cable is in the accessory kit.

❖ To install the serial communication cable

1. Connect one end of the serial communication cable to an available serial port on the host computer.
2. Connect the other end of the cable to the connector found on the back of the Surveyor MS Pump Plus (see [Figure 7](#)).

System Interconnect Cable

The interconnect cable coordinates the timing of the Surveyor Plus modules during an injection sequence.

The Surveyor Plus LC devices are shipped with two versions of the interconnect cable: a 5-connector version (F5049-010) and a 7-connector version (60053-63034).

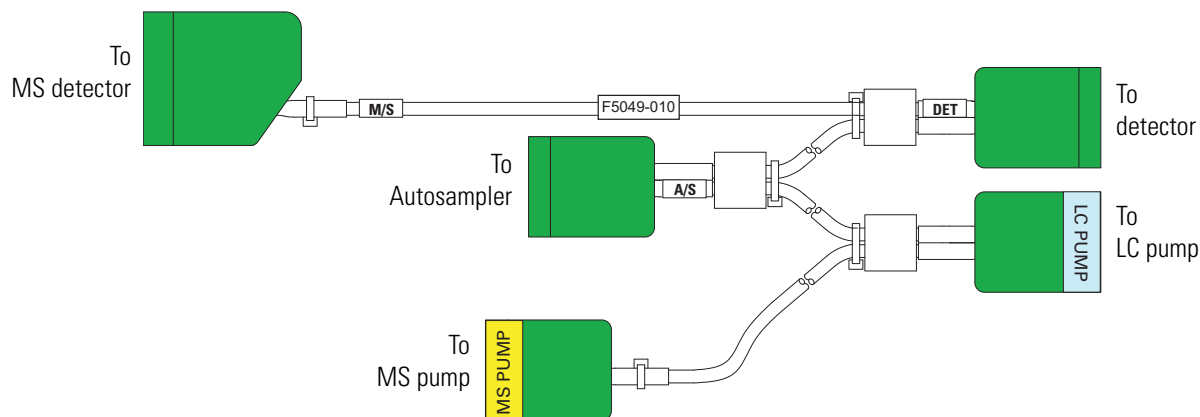
This topic contains the following procedures:

- [Connecting the 5-Connector Interconnect Cable to the LC Devices](#)
- [Connecting the 7-Connector Interconnect Cable to the LC Devices](#)

Connecting the 5-Connector Interconnect Cable to the LC Devices

The 5-combicon connectors of the interconnect cable (P/N F5049-010) are labeled LC PUMP, MS PUMP, A/S, DET, and M/S. See [Figure 8](#).

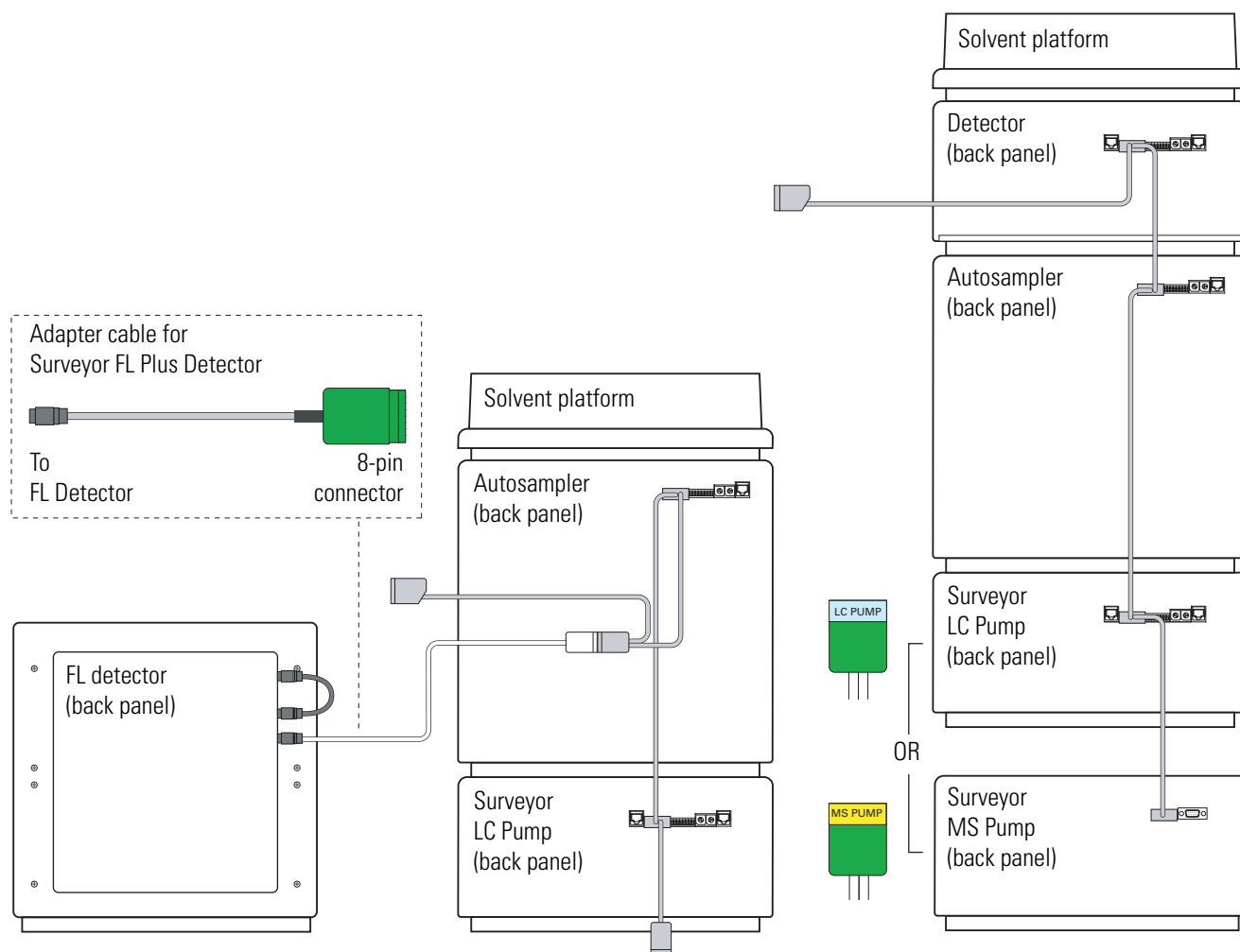
Figure 8. 5-connector interconnect cable, P/N F5049-010



❖ **To connect the LC modules with the 5-connector interconnect cable**

1. To connect the autosampler, plug the A/S connector into the left, 8-pin socket on the back panel of the autosampler. See [Figure 9](#).
2. To connect the MS pump, plug the connector identified by the yellow MS PUMP sticker into the 8-pin receptacle on the back panel of the pump. There is a yellow MS PUMP label above the receptacle.
3. To connect a Surveyor PDA Plus Detector, plug the DET connector into the left, 8-pin socket on the back panel of the detector.
4. To connect a Surveyor FL Plus Detector, plug the DET connector into the adapter cable for the Surveyor FL Plus Detector. Then connect the other end of the adapter cable to the FL detector.

Figure 9. Back panels, showing the connections for the 5-connector interconnect cable



Not drawn to scale

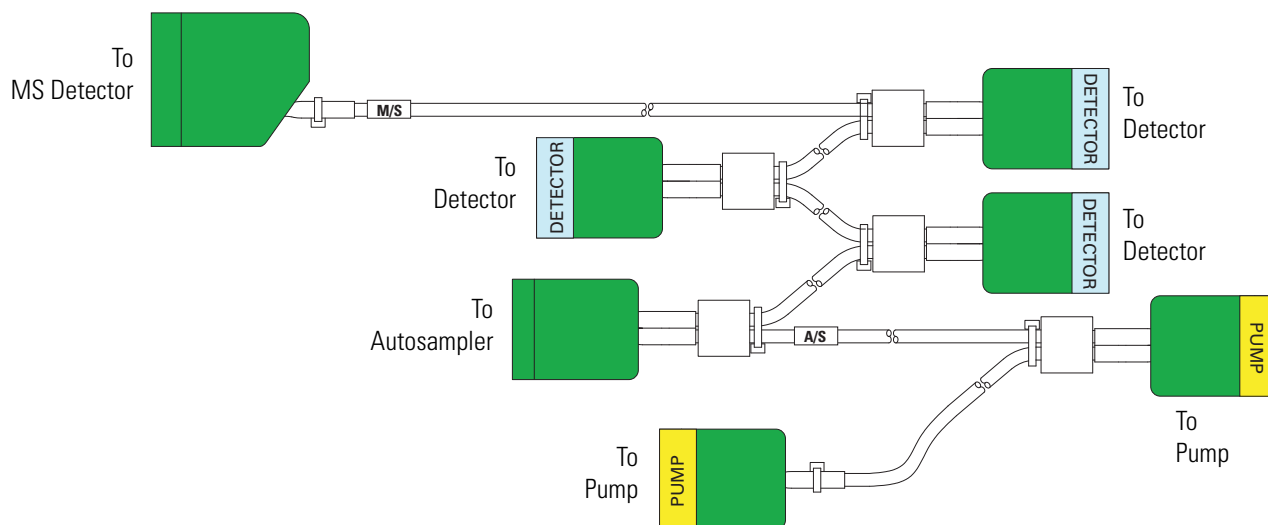
2 Installation

Making the Initial Back Panel Connections

Connecting the 7-Connector Interconnect Cable to the LC Devices

The 7-combicon connectors of the interconnect cable are labeled as follows: three connectors are labeled DETECTOR; two connectors are labeled PUMP; the connector for the A/S is identified by the A/S tag on its adjacent cable; the connector for the MS detector is identified by the M/S tag on its adjacent cable. See [Figure 10](#).

Figure 10. 7-connector version of the system interconnect cable



❖ To connect the LC modules with the 7-connector interconnect cable

1. To connect the Surveyor Autosampler Plus, plug the A/S connector into the left, 8-pin socket on the back panel of the autosampler.
2. To connect a Surveyor PDA Plus Detector, plug the detector connector into the left, 8-pin socket on the back panel of the detector.
3. To connect a Surveyor MS Pump Plus, do the following:
 - a. Connect the end of the adapter cable labeled PUMP to one of the pump connectors of the 7-connector interconnect cable. The adapter cable for the MS pump is shown in [Figure 11](#).
 - b. Plug the end of the adapter cable labeled MS PUMP into the 8-pin socket on the back panel of the MS pump. See [Figure 12](#).

Figure 11. Adapter cable for the Surveyor MS Pump Plus

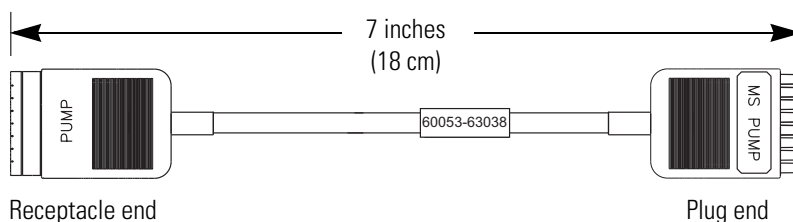
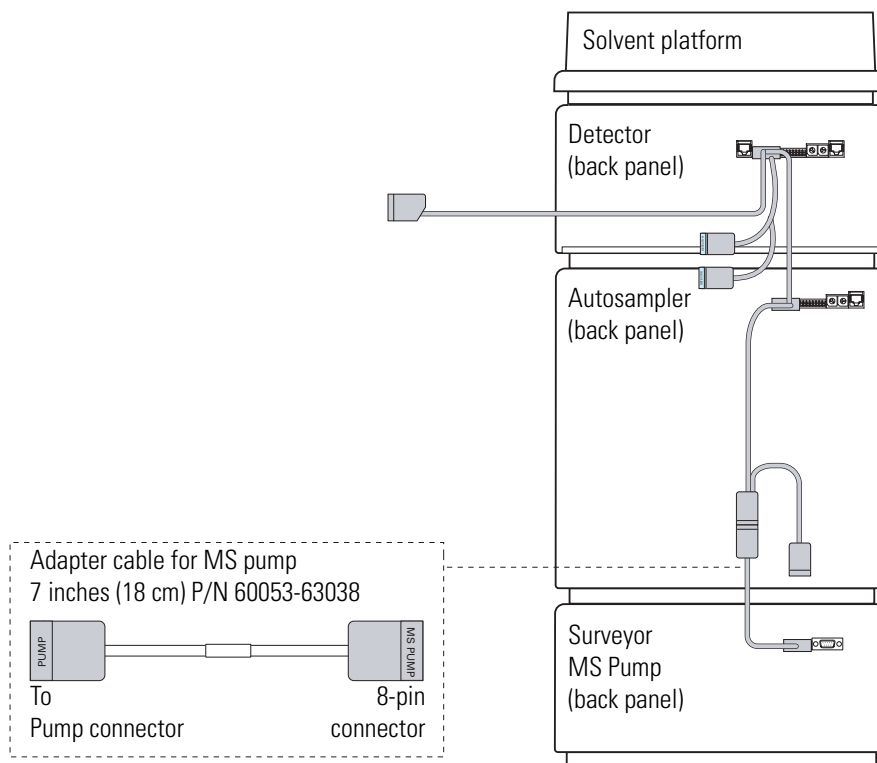


Figure 12. Back panels, showing the connections for the 7-connector interconnect cable

Connecting the Low-Pressure Solvent Lines

You need to connect the low-pressure solvent lines to the Surveyor MS Pump Plus before you operate the pump. There are four low-pressure lines that bring the four solvents to the pump.

Solvents should always be HPLC quality and free of particulate matter. Because the Surveyor MS Pump Plus contains very small, precision-machined parts, it can easily be affected by small particles entering the pump. Aqueous eluents are particularly susceptible to bacterial growth that can cause numerous problems.

To prevent bacteria and/or particulate matter from entering the pump, do the following:

- Filter all solvents thoroughly.
- Discard aqueous mobile phases on a daily basis or add a bactericide such as sodium azide to inhibit bacterial growth (0.65 g/l NaN_3 prevents growth of most kinds of bacteria).



CAUTION Sodium azide is highly toxic and carcinogenic. Take care when disposing of solutions that contain sodium azide because azides can react with heavy metals to form explosive compounds.

2 Installation

Connecting the Low-Pressure Solvent Lines

Air permeates through FEP tubing and can saturate solvents inside the solvent lines that connect the degassing assembly to the solvent proportioning assembly when the system is left static (for example, over a weekend). In this case, purging the lines for a few minutes proves effective.

The Surveyor System Accessory Kit (P/N SRVYR-SYSKT) contains 240 in. of 1/8 in. OD, 1/16 in. ID tubing to cut and use as the solvent inlet lines. Super Flangeless fittings are used to connect these lines to the inlet ports of the vacuum degas assembly of the Surveyor MS Pump Plus. The system accessory kit also contains solvent bottles, solvent bottle caps, and stick on labels for the bottle caps.

To install the solvent inlet lines, perform the following procedures:

- [Step 1: Assembling the Solvent Reservoir Bottles](#)
- [Step 2: Installing the Super Flangeless Fittings](#)

Step 1: Assembling the Solvent Reservoir Bottles

❖ To assemble the solvent reservoir bottles

1. Cut the 1/8 in. OD tubing into four sections of approximately 5-ft. each. Make sure that you cut the ends of the tubing squarely at a 90° angle.
2. Place the label A sticker onto one of the four solvent reservoir bottle caps.
3. Pass one end of the inlet line for solvent A through the solvent reservoir bottle cap labeled A. Terminate the inlet line with one of the solvent reservoir filters included in the system accessory kit. See [Figure 13](#).

Note A Teflon end-of-line filter is supplied with the system accessory kit for this pump. Use only non-metallic filters on the solvent lines.

Figure 13. Solvent reservoir cap assembly



4. Insert the solvent reservoir filter and inlet line into one of the solvent reservoir bottles, and screw the cap onto the solvent reservoir bottle until it is secure.

Note The cap is a two-piece assembly. The upper section (Figure 13) snaps onto a threaded section. The threaded section can be screwed onto the bottle and the upper section can be snapped on after the tubing has been installed, or, if you are replacing existing tubing, the entire cap can be unscrewed from the bottle.

5. Position the bottle in the Surveyor Solvent Platform, allowing the solvent inlet line to hang down along the left side of the Surveyor Plus system.

Step 2: Installing the Super Flangeless Fittings

The Super Flangeless™ fitting consists of three components: a nut, a ferrule, and a stainless steel compression ring (see Figure 14). The compression ring has two sides (see Figure 15). When you place the three components of the fitting on the end of a solvent line, the flattened side of the ring faces the nut and the angled side of the ring faces the ferrule.

Figure 14. Super Flangeless fitting components

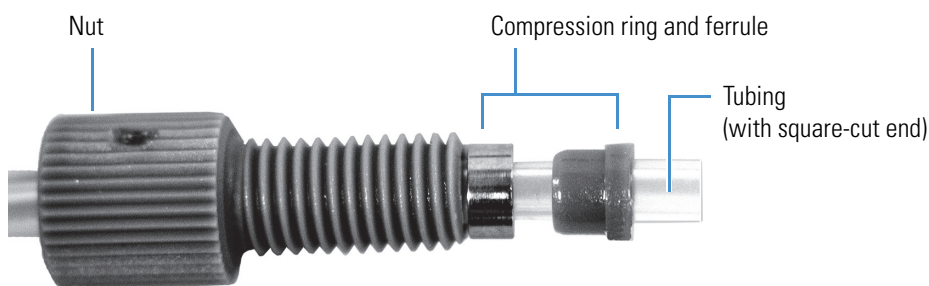
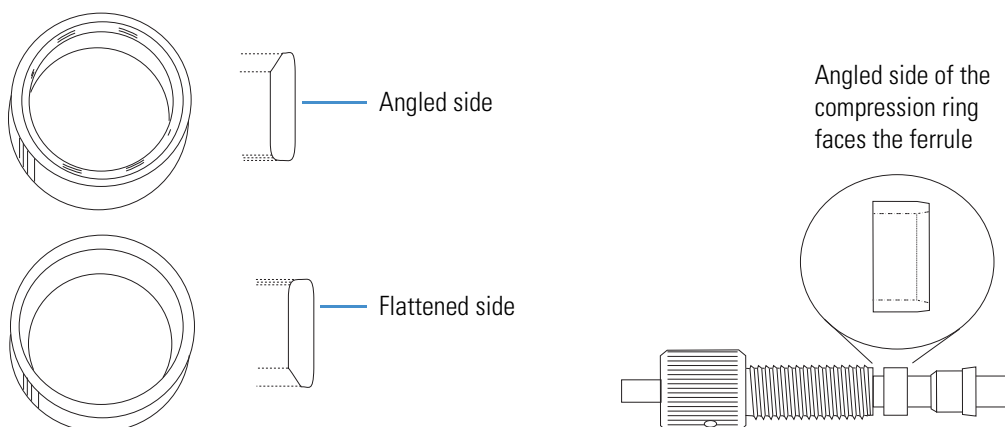


Figure 15. Drawing of the compression ring sides



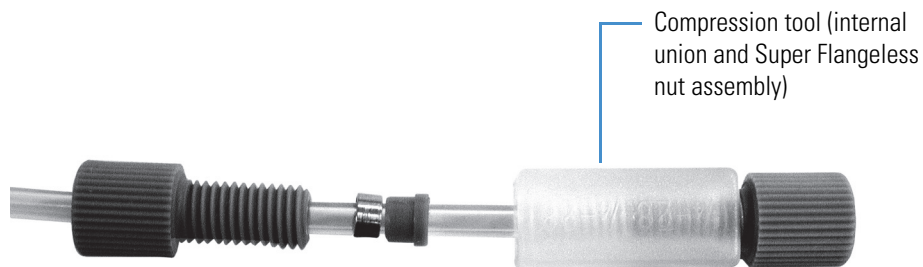
2 Installation

Connecting the Low-Pressure Solvent Lines

❖ To install a Super Flangeless fitting

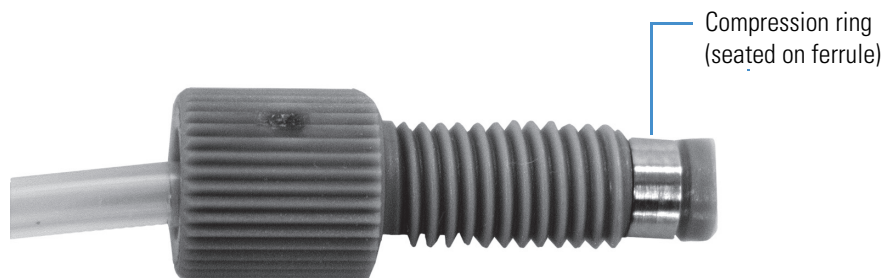
1. Slide the nut onto the end of the tubing.
2. With its flattened side facing toward the nut, slide the compression ring onto the end of the tubing.
3. With its narrow end facing the compression ring, slide the ferrule onto the end of the tubing.
4. Swage the fitting onto the tubing (see [Figure 16](#)):
 - a. Create a compression tool by screwing a Super Flangeless nut into one end of a 1/4 inch × 28 thread internal union.
 - b. Insert the tubing with the fitting assembly into the other end of the union.
 - c. Hold the tubing to the bottom of the tool while tightening down the nut.

Figure 16. Using a compression tool to swage the fitting onto the tubing

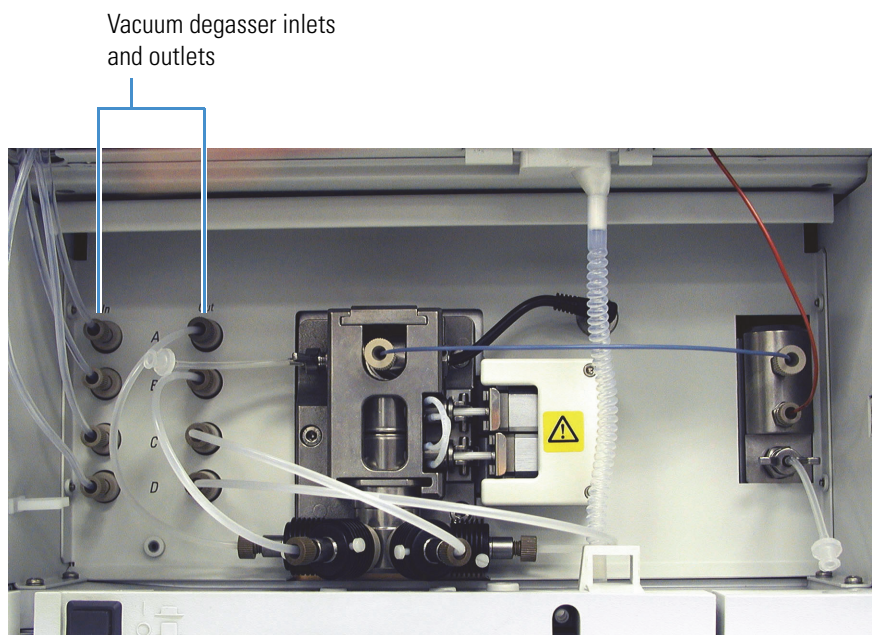


- d. Unscrew the swaged fitting and verify the following (see [Figure 17](#)):
 - The end of the square-cut tubing is flush with the end of the ferrule.
 - The steel compression ring is seated over the ferrule.

Figure 17. View of fitting swaged onto tubing



5. Insert the tubing with the properly swaged fitting into inlet port A, and then tighten the nut fingertight. Be careful not to cross-thread the fitting, which can cause solvent leakage (see [Figure 18](#)).
6. Pass the solvent inlet line through the openings along the left side of each Surveyor Plus module to complete the installation.
7. Repeat the above steps for the B, C, and D solvent inlet lines.

Figure 18. Vacuum degas connection

Connecting the Pump to the Surveyor Autosampler Plus

The solvent outlet tubing is a 1/16 in. stainless steel tube that is used to deliver mobile phase from the Surveyor MS Pump Plus to the Surveyor Autosampler Plus. The tubing, which is pre-installed on the autosampler inlet, has an internal diameter of 0.020 in. A 1/16 in. stainless steel nut and PEEK ferrule, Fingertight fitting is included to connect the tubing to the outlet fitting of the pulse dampener of the MS pump.

❖ To connect the MS pump to the inlet tubing of the autosampler

1. Place the Surveyor Autosampler Plus on top of the Surveyor MS Pump Plus. Locate the autosampler inlet tubing (extending out of the column oven) to be connected to the pump.
2. Carefully bend the Surveyor Autosampler Plus inlet tubing so that it passes through the access port found on the bottom-right side of the Surveyor Autosampler chassis.
3. Connect the tubing to the outlet of the pulse dampener using the 1/16 in. stainless steel nut and PEEK ferrule (P/N 00101-18088) included in the Surveyor MS Pump Plus accessory kit. Be careful not to over-tighten the fitting.

Note To decrease the gradient delay volume of your system, bypass the tubing behind the column oven. This tubing allows the mobile phase to equilibrate to the requested column oven temperature, but it also adds 250 μL of gradient delay volume. To bypass this tubing, disconnect it from port 5 of the injection valve, and then connect the pulse dampener outlet directly to this port.

2 Installation

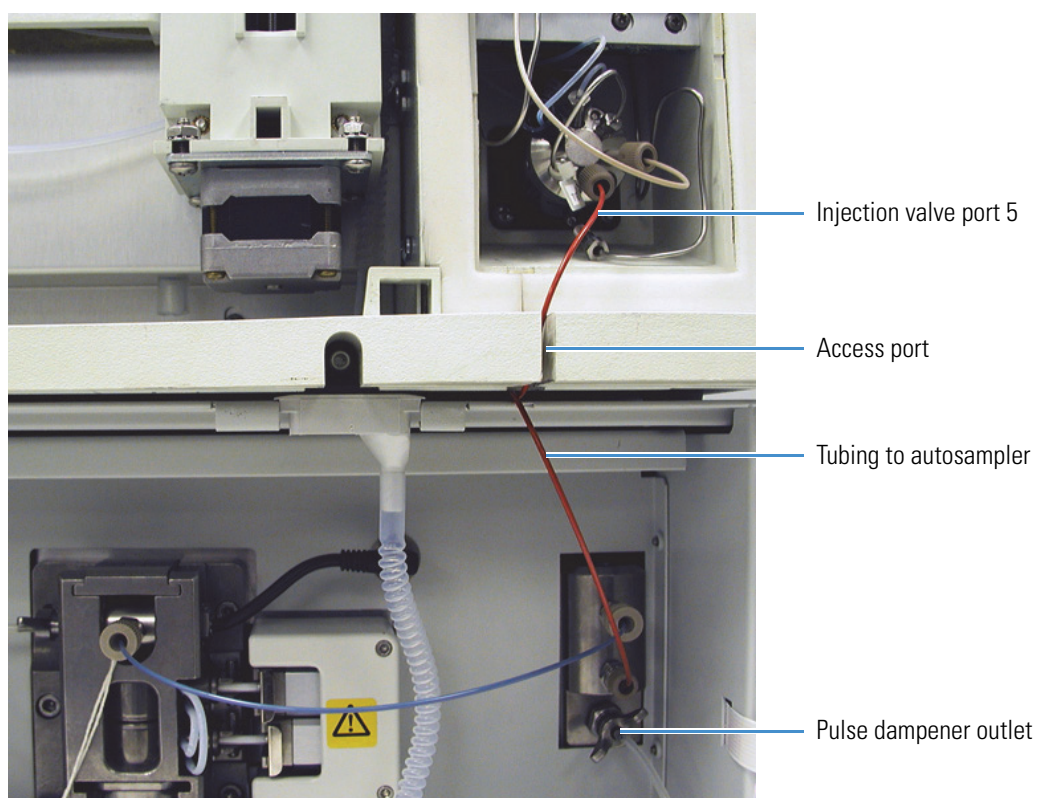
Connecting the Pump to the Surveyor Autosampler Plus

❖ To connect the pump directly to the autosampler's injection valve

1. Place the Surveyor Autosampler Plus on top of the Surveyor MS Pump Plus.
2. Disconnect the stainless steel tubing from port 5 of the Autosampler's injection valve.
3. Use 0.005 in. ID PEEK tubing or the 30 cm (12 in.) l, 0.010 in. ID, stainless steel tubing (P/N A0941-010) provided in the autosampler accessory kit, a Valco injection valve fitting, and the 1/16 in. stainless steel nut and PEEK ferrule (P/N 00101-18088) included in the Surveyor MS Pump Plus accessory kit to connect the outlet of the pulse dampener directly to port 5 of the injection valve. See [Figure 19](#).

The Surveyor MS Pump Plus is now properly connected.

Figure 19. Solvent outlet tubing connection to autosampler



Powering On for the First Time

After the solvent lines and tubing are installed, the pump is ready to be powered on.

❖ To turn on the pump

1. Ensure that the power switch at the front of the unit is in the Off position (released or Out-position).
2. Attach the power cord to the power entry module on the back panel of the pump and connect to the power source (Figure 7).
3. Push the power On button to engage it (In-position).
4. Note that the power indicator turns solid green (see Figure 6 on page 6). If it does turn green, see “Surveyor MS Pump Troubleshooting Guide” on page 68. Call your Thermo Fisher Scientific service representative if you require assistance.

Configuring the Data System PC COM Port

Communication with the Surveyor MS Pump is established through an RS232 connection.

❖ To configure the COM port



1. From the Windows taskbar, choose **Start > Control Panel**.
2. Double-click the **System** icon to open the System Properties dialog box.
3. Click the **Hardware** tab, and then click **Device Manager**.
4. Double-click **Ports (COM & LPT)**. The available ports are displayed below Ports (COM & LPT) in the Device Manager list.
5. Double-click **Communication Port (COM1)** to display the Communication Port (COM1) Properties dialog box.
6. Click the **Port Setting** tab.
7. Set the configuration parameters
 - Bits per Second: 19200
 - Data Bits: 8
 - Parity: none
 - Stop Bits: 1
 - Flow Control: none
8. Click **OK** to save the changes and close the Communication Port (COM1) Properties dialog box.

9. Close the Device Manager window.
10. Click **OK** to close the System Properties dialog box.
11. Restart the computer to enable the new settings.

Preparing the Pump for Operation

After you connect the solvent lines to their respective degas inlets, you are ready to prepare the pump for operation. To prepare the pump for operation, first configure the MS pump to communicate with the Xcalibur data system. Then purge the air out of the solvent lines, and prime the pulse dampener.

This section contains the following topics:

- [Specifying the Instrument Configuration Options](#)
- [Purging Air from the Solvent Lines](#)
- [Priming the Pulse Dampener](#)

Specifying the Instrument Configuration Options

You can use the Surveyor MS Pump Plus (or Surveyor MS Pump) with the Surveyor Autosampler Plus or the CTC PAL autosampler (Thermo PAL driver). You can also set up a system consisting of two Surveyor pumps.

This topic describes the Xcalibur instrument configuration options for the Surveyor MS Pump Plus as well as the configuration options for the Surveyor Autosampler Plus that are required to ensure that the autosampler recognizes the output signals from the pump.

Note The user interface for the Instrument Configuration application compatible with Xcalibur 2.1 or higher has a slightly different appearance than the Instrument Configuration application compatible with earlier Xcalibur versions.

To specify the instrument configuration options, follow these procedures:

- [Opening the Instrument Configuration Application](#)
- [Adding the Surveyor MS Pump or Surveyor MS Pump Plus to the Instrument Configuration](#)
- [Configuring the Signal Polarity for the Surveyor AS](#)

Opening the Instrument Configuration Application

Use the Instrument Configuration (Thermo Foundation Instrument Configuration) window to specify the system hardware components.

❖ To open the Instrument Configuration window for systems using Xcalibur 2.1

From the Windows taskbar, choose **Start > All Programs > Thermo Foundation > Instrument Configuration**.

The Thermo Foundation Instrument Configuration window appears.

❖ To open the Instrument Configuration window for systems using Xcalibur 2.0.x or lower

- From the Windows taskbar, choose **Start > All Programs > Xcalibur > Instrument Configuration**.
- Or, double-click the Xcalibur Instrument Configuration icon on the Windows desktop.

The Instrument Configuration window appears.

Adding the Surveyor MS Pump or Surveyor MS Pump Plus to the Instrument Configuration

❖ To add the pump to the Instrument configuration

1. Open the Instrument Configuration window.
2. In the Available Devices list, double-click the icon for the pump.

A copy of the icon appears in the Configured Devices list.

3. In the Configured Devices list, double-click the pump icon.

The Surveyor MS Pump Plus (or Surveyor MS Pump) dialog box appears.

Configuring the Signal Polarity for the Surveyor AS

During an injection sequence, the Surveyor MS Pump Plus issues a Pump Ready signal and an Injection Hold Release signal to the autosampler.

When you configure the Surveyor Autosampler Plus, select the Pump ready active high check box and the Injection hold release active high check box. If you do not select these two check boxes, the autosampler cannot recognize the input signals from the pump and does not begin an injection sequence.

After you submit a run, the Surveyor Autosampler Plus waits for a Pump Ready signal from the pump. The pump issues this signal after its pressure transducer monitors a stable backpressure. When the autosampler receives this signal, it begins an injection sequence.

2 Installation

Preparing the Pump for Operation

After the autosampler begins the injection sequence, it waits for an Injection Hold Release signal from the pump. The pump issues this release signal when its piston cam returns to the start position. When the autosampler receives the Injection Hold Release signal, it switches the injection valve to the inject position, allowing the contents of the sample loop to be swept into the mobile phase stream from the pump.

Note In Xcalibur, the instrument setup method for the pump contains a custom stability limit parameter. This parameter allows the user to define the acceptable level of pressure stability that must be attained before the pump issues the Pump Ready Signal to the autosampler.

❖ To configure the input signal polarities for the Surveyor Autosampler

1. If it is not already open, open the Instrument Configuration window (see “Opening the Instrument Configuration Application” on page 23).

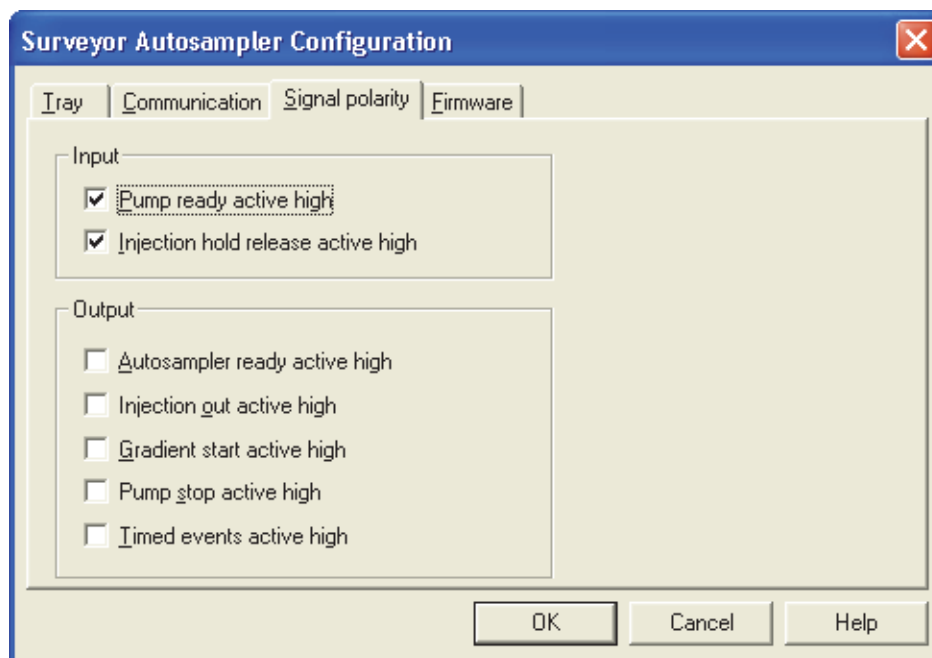
2. In the Configured Devices list, double-click the **Surveyor AS** icon.

The Surveyor Autosampler dialog box appears.

3. Click the **Signal polarity** tab.

The Signal polarity page appears (see Figure 20).

Figure 20. Surveyor Autosampler Configuration dialog box



4. In the Input area, make sure that the **Pump ready active high** and **Injection hold release active high** check boxes are selected.

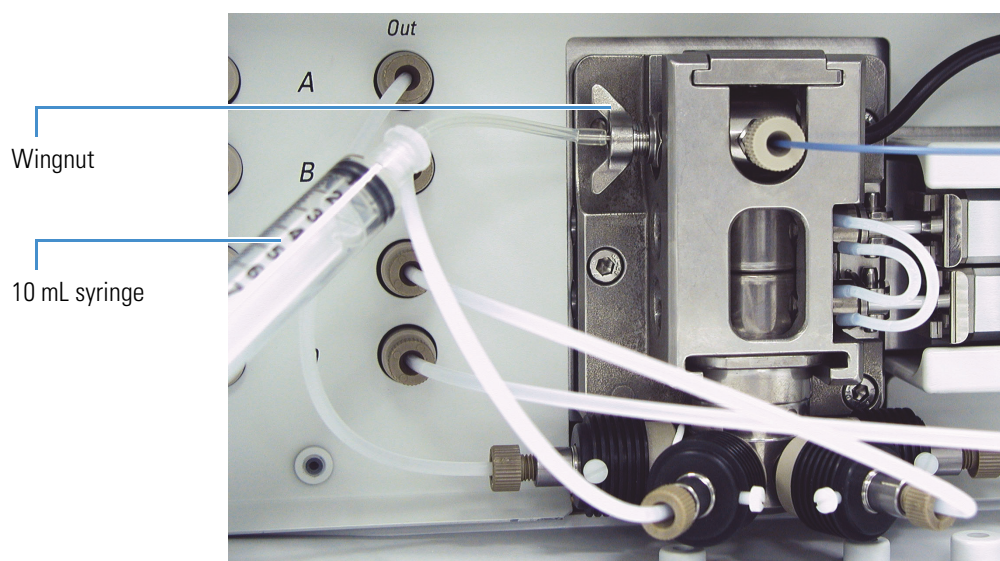
Purging Air from the Solvent Lines

After you fill the solvent bottles, connect the solvent lines to the degasser, and configure the MS pump to communicate with Xcalibur, prepare the pump for operation by removing the air from the solvent lines.

❖ To remove air from the solvent lines

1. Insert the tip of the 10 mL syringe into the tubing that is connected to the wingnut on the left side of the LDA. See [Figure 21](#).
2. Open the wingnut by turning it counterclockwise.

Figure 21. Purging the solvent lines



3. Turn on the pump power.
4. Open the Direct Control dialog box for the pump:
 - a. Open Xcalibur.
 - b. On the Roadmap view, click the **Instrument Setup** button. Or, from the menu bar, choose **GoTo > Instrument Setup**.
 - c. In the view bar, click the **Surveyor MS Pump Plus** (or **Surveyor MS Pump**) button.
 - d. Choose **Surveyor MS Pump Plus** (or **Surveyor MS Pump**) > **Direct Control** from the Instrument Setup window to open the Direct Control dialog box.
5. In the Direct Control dialog box (see [Figure 22](#)), do the following:
 - a. Select the **Take pump under control** check box.
 - b. Type percentages in the Inlet boxes for the solvent lines through which you want to draw solvent.
 - c. In the Flow box, type **2000**.



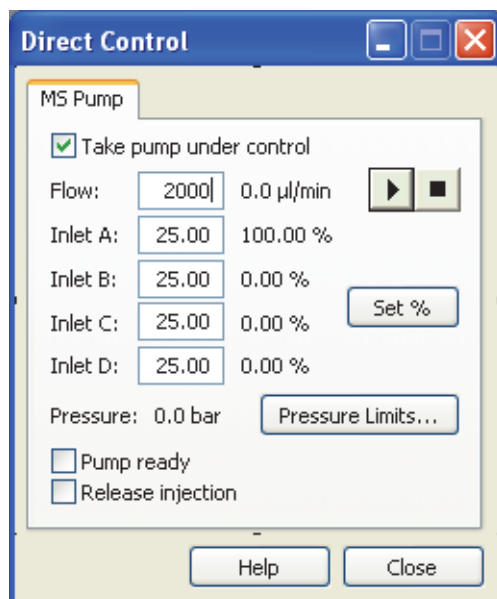
2 Installation


Preparing the Pump for Operation

- d. Click the **Start Run** button  to start the pump flow.

Because each 1.5 m (5-ft.) section of tubing holds approximately 3 mL of solvent, you need to stop the pump flow and empty the syringe periodically.

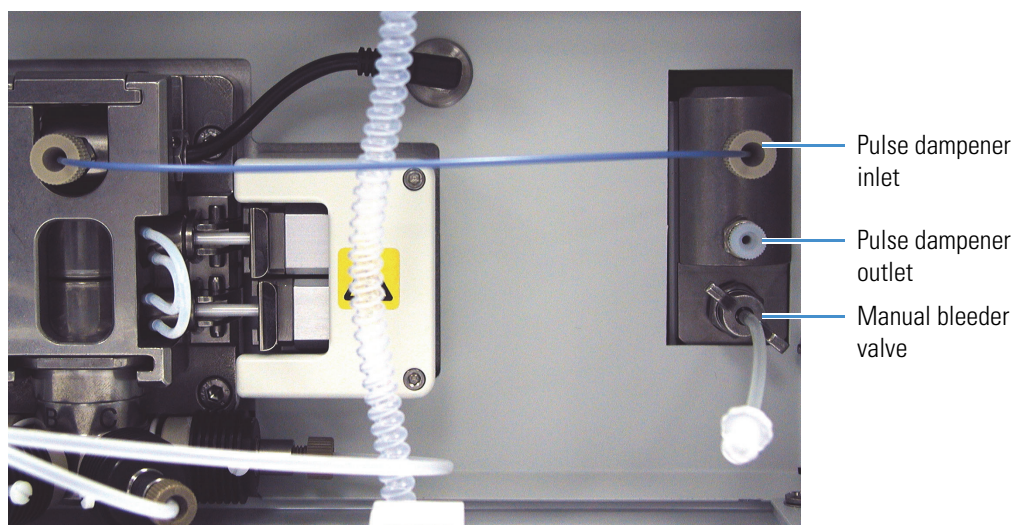
Figure 22. Direct Control dialog box for the pump



6. After you finish purging the solvent lines, click the **Stop** button  to stop the pump flow.
7. Turn the LDA wingnut clockwise to close the venting valve.

Priming the Pulse Dampener

Priming the pulse dampener involves filling the dampening loop of the pulse dampener with a solvent such as methanol or isopropanol. After the loop is filled with solvent, it can effectively dampen pressure pulsations from the rest of your system. Closing the manual bleeder valve after you fill the loop shuts the loop off from the flow path. For standard operation, reprime the pump when you notice an increase in the pump pulsation. If you accidentally open the manual bleeder valve on the front of the pulse dampener while you are pumping an acidic or a buffered mobile phase, immediately flush the buffer out of the loop, and then prime the pulse dampener. See [Figure 23](#).

Figure 23. Pulse dampener assembly

Note Do not fill the pulse dampener with an aggressive acid or a buffered solution. The recommended filling solvents are methanol, acetonitrile, or isopropyl alcohol. If you accidentally fill the pulse dampener with an acidic or buffered solvent, flush the loop with a miscible solvent, and then refill it with methanol, acetonitrile, or isopropyl alcohol.

❖ **To prime the pulse dampener**

1. Fill a solvent reservoir with particulate-free, HPLC grade methanol.
2. Connect an HPLC column or a flow restrictor to the pulse dampener outlet.
3. Insert the tip of the 10 mL syringe into the tubing that is connected to the manual bleeder valve of the pulse dampener, and then open the valve by turning it counterclockwise.
4. Set the flow rate to 1 mL/min.
5. Fill the loop completely, to expel any air that might be trapped in the dampener loop.
6. Set the pump flow to a rate that is appropriate for your system.
7. Close the manual bleeder valve of the pulse dampener by turning the valve clockwise.

The pulse dampener loop is now filled. Unless you need to refill the dampener loop of the pulse dampener with fresh solvent, keep the manual bleeder valve closed.

Diagnostics

This chapter contains the following diagnostic procedures that can be performed on the Surveyor MS Pump Plus.

Contents

- [Monitoring Pump Pulsation](#)
- [Testing the Pump Proportioning](#)
- [Determining the Delay Volume of the Pump](#)

Monitoring Pump Pulsation

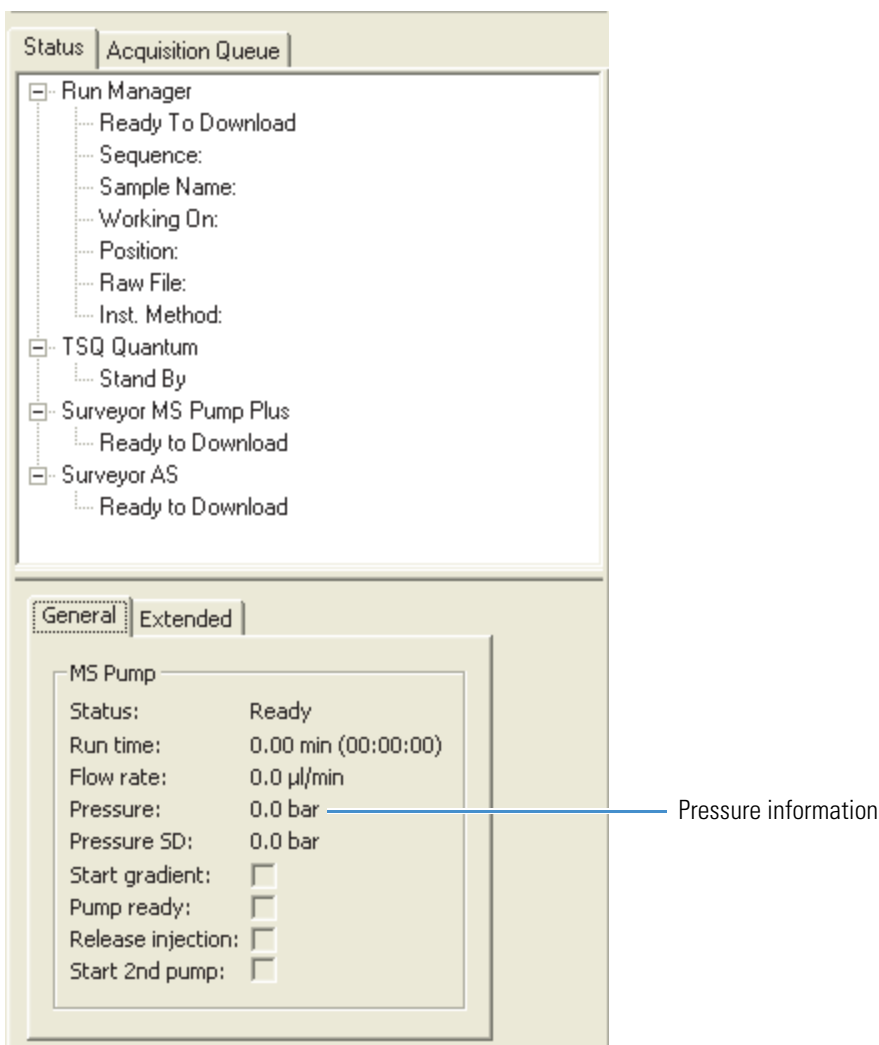
The pump pulsation should be less than 1% at backpressures greater than 100 bar (1450 psi).

Note The Surveyor MS Pump Plus pressure fluctuation often appears to be higher than that for conventional pumps, because the Surveyor MS Pump Plus does not use electronic pulse damping.

The pressure pulsation can be directly monitored through the Xcalibur data system. You can find the pressure information in the Status page of the Information view as shown in [Figure 24](#). This view is normally displayed on the left side of the Home Page window.

If the Info view is not displayed, it has been turned off. From the Home Page window, choose **View > Info View** to toggle the Information view On and Off. Then, click the Status tab to display the Status page.

Figure 24. Status page of the Information view



Trapped air bubbles, contaminated proportioning valves, or contaminated check valves can cause excessive pump pulsation. If the pulsation of the pump is greater than 1% at pressures greater than 100 bar, purge air out of the piston chambers as described on [page 31](#), and clean the proportioning valves and the check valves as described in “[Cleaning the LDA Components without Disassembling the LDA](#)” on [page 38](#).

If you cannot reduce the variation in backpressure by performing these procedures, disassemble and clean the LDA and replace the primary piston seals as described in “[Replacing the Primary Piston Seals](#)” on [page 49](#).

Testing the Pump Proportioning

You need to test the proportioning accuracy of the pump periodically by running a step-gradient.

Two eluents, methanol and spiked methanol, are used in this procedure. The absorbance reading for the 100% spiked methanol solution should be in the 600 to 800 mAU range.

- If your detector contains a standard 10 mm flow cell, prepare the spiked methanol solution by adding 7 mg of ethyl paraben (P/N A4904-010) to 1 L of HPLC grade methanol.
- If your detector contains a 50 mm LightPipe flow cell, prepare the spiked methanol solution by adding 1.4 mg of ethyl paraben to 1 L of HPLC grade methanol.

The step gradient program is shown in [Table 2](#). Use the following conditions to run the step gradient:

- Flow: 200 μ L/min
- Pressure: approximately 100 bar by restriction
- Eluents: A: methanol
B: spiked methanol
- Sub-proportioning: Off
- Steps: 10 min hold, 20% increments
- Detection: UV (254 nm)

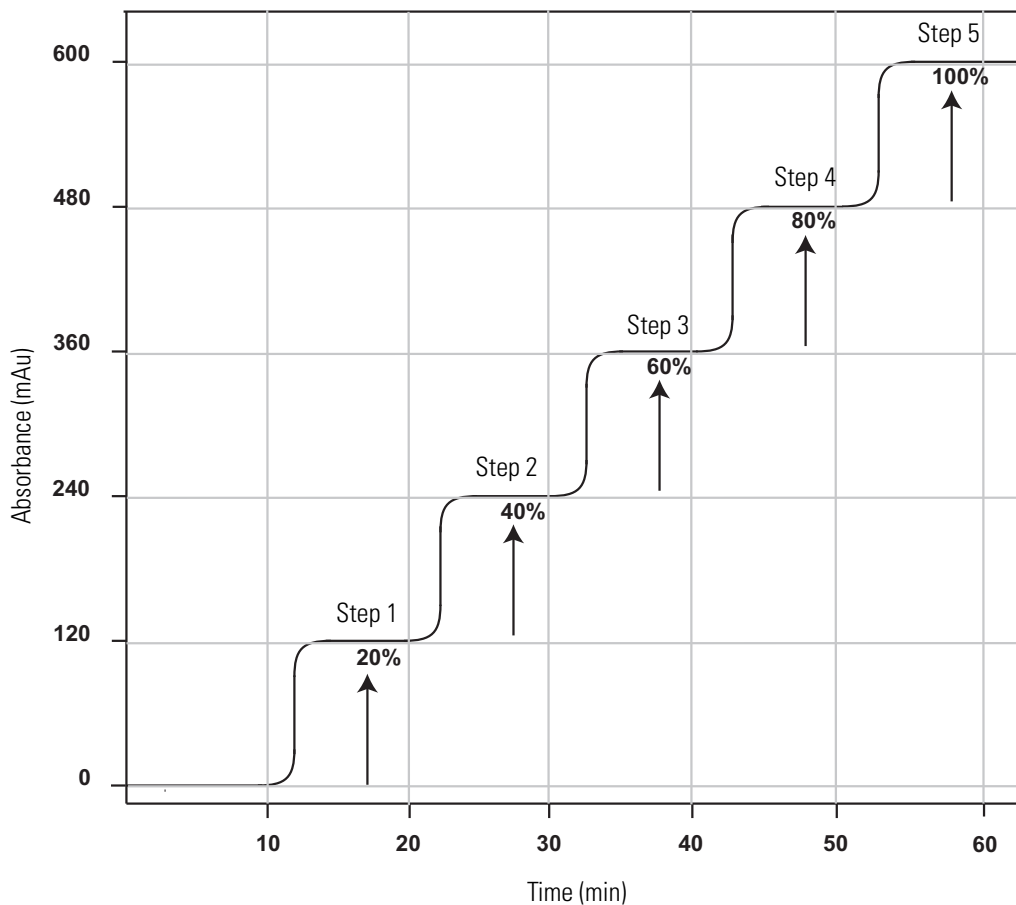
Table 2. Gradient Program

Time (min)	%A	%B
0.00	100.0	0.0
10.0	100.0	0.0
10.01	80.0	20.0
20.0	80.0	20.0
20.01	60.0	40.0
30.0	60.0	40.0
30.01	40.0	60.0
40.0	40.0	60.0
40.01	20.0	80.0
50.0	20.0	80.0
50.01	0.0	100.0
60.0	0.0	100.0

Plumb your system to bypass the autosampler. Use the tubing supplied with the Surveyor system (P/N 803260, Inlet tubing, with insulation, PEEK 1/16 in. OD × .005 in. ID, Red) to connect the pump directly to the detector. Use a flow restrictor to maintain a backpressure greater than 40 bar.

The resulting gradient profile should contain five steps of equal height. See [Figure 25](#). Inaccuracies in the step heights indicate proportioning errors.

Figure 25. Step gradient profile for testing compositional accuracy



Step heights should be within $\pm 1\%$ of the theoretical values. In the step gradient shown in [Figure 25](#), the absorbance value for 100% spiked solvent is 600 mAU (600 mAU – 0 mAU). Therefore, the theoretical values for the 20%, 40%, 60%, and 80% steps are 120, 240, 360, and 480 mAU, respectively. A deviation greater than 6 mAU for any of the steps would be a failing result. If the pump passes both the pulsation and proportioning tests, it is working properly.

If the pump fails the pulsation test or the proportioning test, try to restore proper performance by cleaning the proportioning valves as described in [“Cleaning the LDA Components without Disassembling the LDA”](#) on [page 38](#). If this fails to restore proper pump performance, disassemble and clean the LDA and replace the primary seals as described in [“Replacing the Primary Piston Seals”](#) on [page 49](#).


Determining the Delay Volume of the Pump

The gradient delay volume of your pump is an important variable if you are developing gradient methods for use on multiple LC systems.

The gradient delay volume of your Surveyor MS Pump Plus consists of the wetted components that lie in the flow path between the solvent proportioning valves of the pump and the outlet of the pulse dampener.

Note The tubing located behind the built-in column oven of the Surveyor Autosampler contains approximately 250 μL of gradient delay volume.

❖ To determine the gradient delay volume of your MS pump

1. Set up your system:
 - a. Bypass the autosampler by connecting the pump directly to the detector with tubing of known volume. Each inch of the inlet tubing to the Surveyor PDA detector (P/N 803260, Inlet tubing, with insulation, PEEK, 1/16 in. OD \times 0.005 in. ID, Red) contains 0.32 μL of gradient delay volume.
 - b. Prepare a spiked methanol solution as specified in [“Testing the Pump Proportioning”](#) on [page 31](#).
 - c. Fill one solvent reservoir bottle with 100% methanol.
 - d. Fill one solvent reservoir bottle with the spiked methanol solution.
 - e. Purge the solvent lines until the old solvents are replaced with the new solvents.
2. Create an instrument method:
 - a. In the General page of the Surveyor MS Pump Plus Instrument Setup view, type the solvent names.
 - b. In the Gradient page of the Surveyor MS Pump Plus Instrument Setup view, add a time line for 0.01 minutes. Switch the mobile phase solvent to the spiked methanol solvent at this time point. Enter a flow rate of 200 $\mu\text{L}/\text{min}$ for both the 0.00 min time line and the 0.01 min time line.
 - c. Program your detector to collect one channel at 254 nm for a period of 5 minutes.
 - d. Save the method.
3. Equilibrate the system until the baseline stabilizes:
 - a. In the Direct Control dialog box, take the pump under control, and set the parameters to 100% methanol and a 200 $\mu\text{L}/\text{min}$ flow rate.
 - b. Click the Start button .
4. Create a one-line sequence containing the acquisition method.
5. Run the sequence.

3 Diagnostics

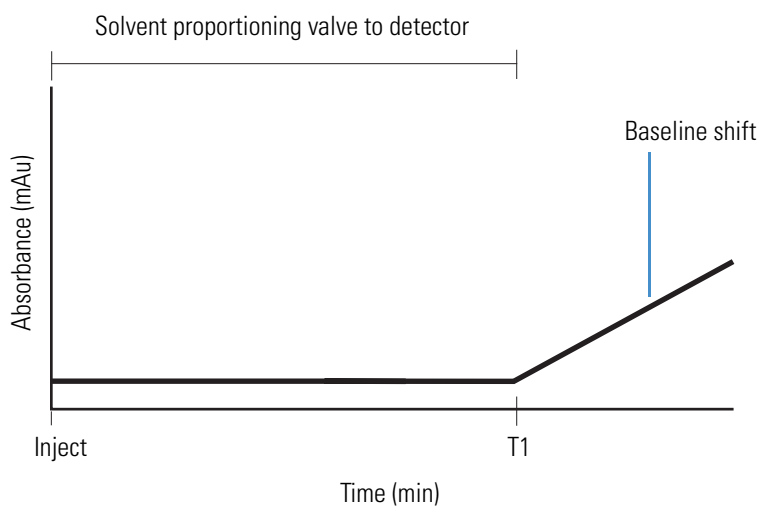
Determining the Delay Volume of the Pump

6. Calculate the gradient delay volume of the pump:
 - a. Determine the total gradient delay volume by multiplying the resulting gradient delay time by the flow rate.
 - b. Determine the gradient delay volume of the pump by subtracting the volume of the tubing connecting the pump to the detector from the total gradient delay volume.

The results of your run should look like the chromatogram shown in [Figure 26](#). If the start of the baseline shift occurs at 0.45 minutes, then the gradient delay volume of the pump is calculated as follows:

$$\begin{aligned}\text{Gradient Delay Volume of Pump} &= (T1 \times \text{Flow Rate}) - \text{tubing volume} \\ &= (0.45 \text{ min} \times 200 \text{ } \mu\text{L}/\text{min}) - 8 \text{ } \mu\text{L} \\ &= 82 \text{ } \mu\text{L}\end{aligned}$$

Figure 26. Graphical display of the baseline shift that occurs as the spiked mobile phase reaches the detector



Maintenance

To properly maintain your pump, monitor its performance. Routinely examine the pump for leaks, and monitor the variation in pump pulsation. In addition, periodically test the proportioning accuracy of the pump. To lengthen the lifespan of the seals, always flush the piston guide bushings with distilled water after pumping buffered solutions.

If the pump pulsation is too high, try removing air from the LDA and cleaning the LDA components. If you detect a leak, disassemble the LDA, clean its components, and replace the primary piston seals.

Contents

- [Flushing the Piston Guide Bushings](#)
- [Purging Air from the Piston Chambers](#)
- [Cleaning the LDA Components without Disassembling the LDA](#)
- [Servicing the LDA](#)
 - [Removing the LDA from the Pump](#)
 - [Disassembling the LDA](#)
 - [Cleaning the LDA Components](#)
 - [Replacing the Primary Piston Seals](#)
 - [Reassembling the LDA](#)
 - [Reinstalling the LDA](#)
- [Servicing the Proportioning Valves](#)
- [Replacing the Solvent Reservoir Filter](#)

Flushing the Piston Guide Bushings

As the pistons move back and forth within the piston chambers, a small quantity of mobile phase leaks behind the primary piston seals. With buffered mobile phases, a precipitate (also known as creep or salt build-up) forms as the liquid fraction evaporates. The abrasion caused by this precipitate can scratch the piston and shorten the life of the piston seals.

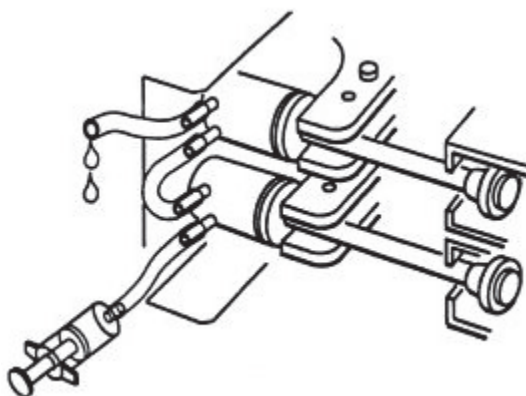
To lengthen the lifespan of the primary piston seals and the pistons, always rinse the piston guide bushings with distilled water after pumping buffered solutions. See [Figure 27](#).

The piston guide bushings are supplied with two pieces of standard 1/8 in. × 1/16 in. PTFE tubing.

❖ To flush the piston guide bushings

1. If the outer piece of PTFE tubing has not been cut, cut it through the middle.
2. Connect a syringe filled with distilled water to the bottom piece of outer tubing and gently flush the water through the system.

Figure 27. Flushing the piston guide bushings



Purging Air from the Piston Chambers

Air trapped in the piston chambers causes an excessive variation in the backpressure.

❖ To purge air out of the piston chambers

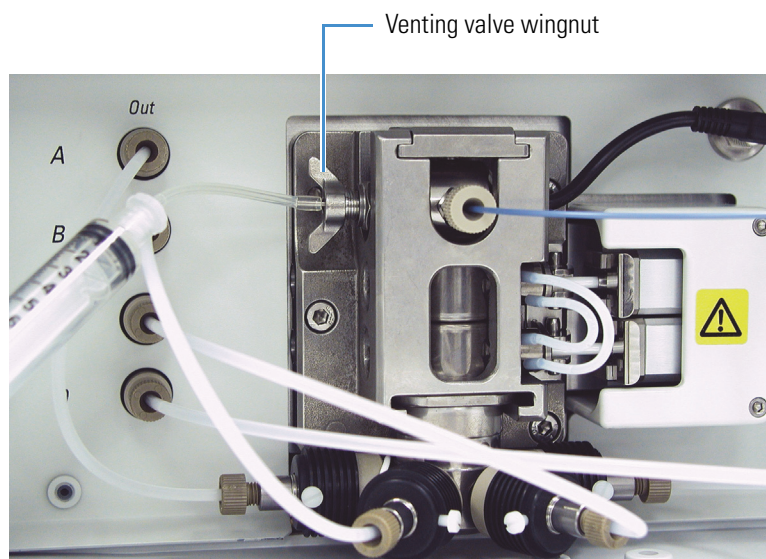
1. Turn off the solvent flow from the pump.



CAUTION Opening the venting valve when the pump is in operation can damage the pressure transducer. Before you open the venting valve, turn off the solvent flow and wait for the system backpressure to dissipate.

2. Open the venting valve by turning the wingnut on the left side of the LDA counterclockwise (see [Figure 28](#)).

Figure 28. Purging air from the piston chambers



3. Insert the 10 mL syringe, which is supplied in the accessory kit, into the tubing attached to the venting valve, and then pump 100% methanol or 100% acetonitrile through the LDA at a flow rate of 200 $\mu\text{L}/\text{min}$ for a period of 10 to 15 minutes.
4. Close the venting valve and return to your initial conditions.
5. Check the pressure variation to determine if performing this procedure remedied the pressure stability problem.

Note Purging the pump at a low flow rate with 100% organic solvent works much better than purging the pump at a high flow rate with (partly) aqueous eluents.

Note All HPLC pumps require proper eluent degassing. The Surveyor MS Pump Plus has a built-in vacuum degasser. Plumbing the pump to bypass this degasser introduces a potentially unacceptable level of air into the piston chambers.

Cleaning the LDA Components without Disassembling the LDA

Contaminated check valves and contaminated proportioning valves can cause an excessive variation in backpressure. To remove the contamination, open the venting valve to bypass the LC column and pump the appropriate flush solvents through the LDA.

❖ To clean the check valves and the proportioning valves

1. Open the venting valve by turning the wingnut on the left side of the LDA counterclockwise (see [Figure 28](#)).

Note Opening the venting valve redirects the mobile phase out the venting valve on the left side of the LDA rather than into the crossover tubing. If you forget to open the venting valve, the flush solvents pass through the pulse dampener and onto your LC column.

2. Set up the system to pump the same solvent on all four channels by connecting all four solvent reservoir lines to one solvent bottle.
3. Attach a 10 mL disposable syringe to the tubing attached to the vent valve. The syringe is supplied in the MS Pump Accessory Kit (see “[MS Pump Accessory Kit](#)” on [page 78](#)).
4. If you were pumping buffered mobile phase through the system, pump HPLC-grade water on all four channels at a flow rate of 200 $\mu\text{L}/\text{min}$ for 30 minutes. Empty the contents of the syringe into a suitable waste receptacle, and then pump an additional 6 mL of water through the LDA.
5. Sequentially pump solvents of increasing polarity through the LDA. The last flush solvent should be miscible with the mobile phase that you plan to use.

Table 3. Polarity indexes of various HPLC solvents

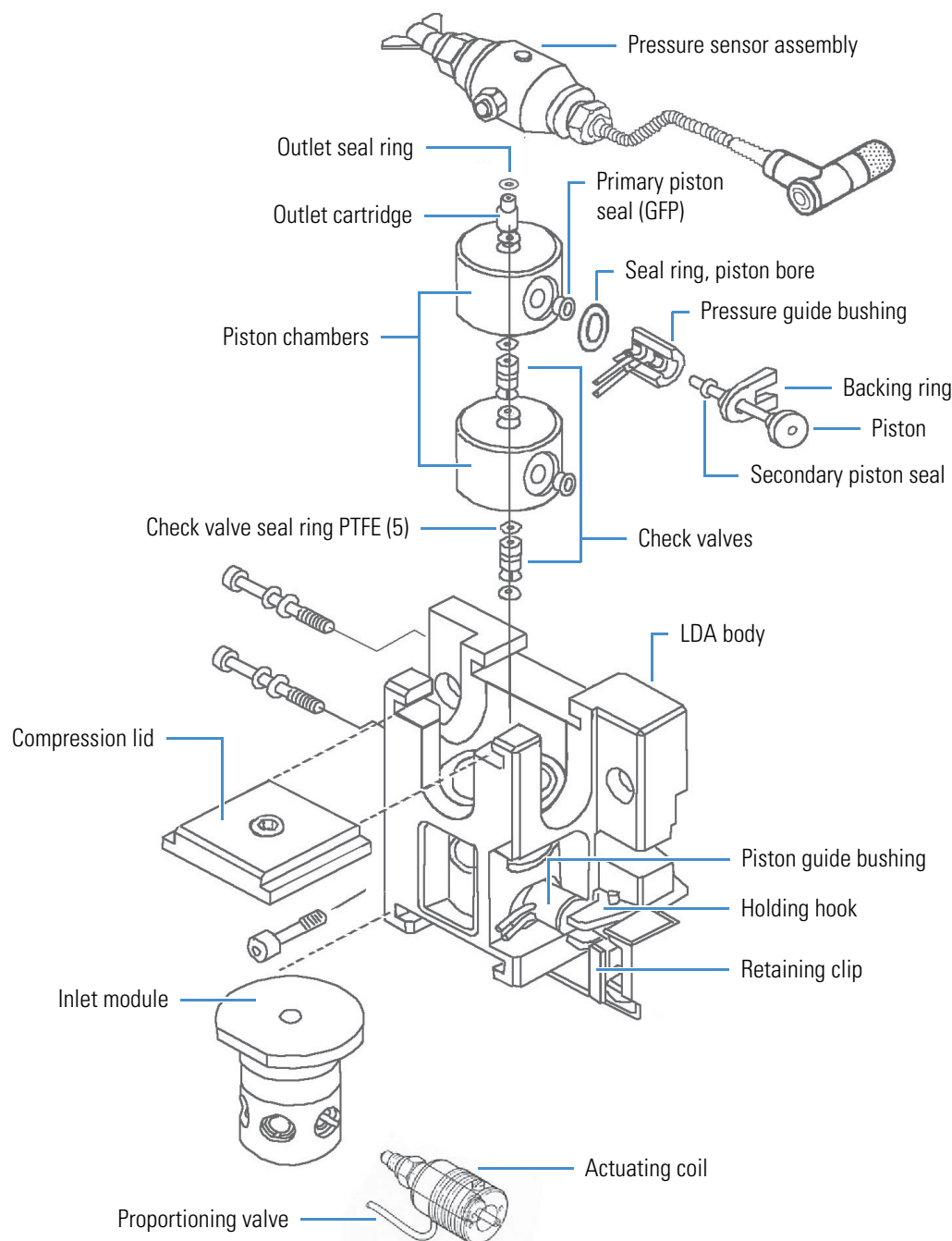
Solvent	Polarity index
Isopropanol	3.9
Methanol	5.1
Acetonitrile	5.8
Water	9.0

6. Close the vent valve.
7. Retest the pressure variation by returning to the initial conditions that displayed the problem. Make sure that the backpressure at these conditions is at least 40 bar.

Servicing the LDA

The liquid displacement assembly (LDA) contains the parts of the pump that are subject to routine wear and tear. These parts are the primary piston seals, the pistons, and the check valves. Occasionally, the secondary piston seals, the check valve seal rings in the top and bottom of the piston chambers, and the outlet seal ring might also require replacement.

Figure 29. Components of the liquid displacement assembly (LDA)



To service the liquid displacement assembly (LDA), follow these procedures:

1. [Removing the LDA from the Pump](#)
2. [Disassembling the LDA](#)
3. [Cleaning the LDA Components](#)
4. [Replacing the Primary Piston Seals](#)
5. [Reassembling the LDA](#)
6. [Reinstalling the LDA](#)

Removing the LDA from the Pump

The must have the following tools to remove the LDA from the pump:

- 4 mm balldriver (with the red screwdriver style handle) (P/N 00725-00032)
- 4 mm Allen wrench (P/N 000725-00034)
- Flat-blade screwdriver

❖ To remove the LDA from the pump

1. Turn off the pump.
2. Remove the crossover tube from the LDA by unscrewing the FingerTight fitting. See [Figure 30](#).
3. Disconnect the pressure sensor cable ([Figure 30](#)) by unscrewing it counterclockwise and pulling it forward.
4. Use the Allen wrench to pry off the retaining clips that secure the pistons to the motor drive arms. Place one end of the wrench behind the lip of the clip, and then push the other end of the wrench toward the pump face. See [Figure 31](#).
5. Raise the solvent filter in each bottle above the solvent level so that only the fluid in the lines drains out. Use a beaker to collect the mobile phase contained in the lines.
6. Unscrew the solvent lines (1/8 in. ID Teflon tubing) connected to the four solvent proportioning valves.

Figure 30. Front of MS pump, showing the crossover tube and the pressure sensor cable

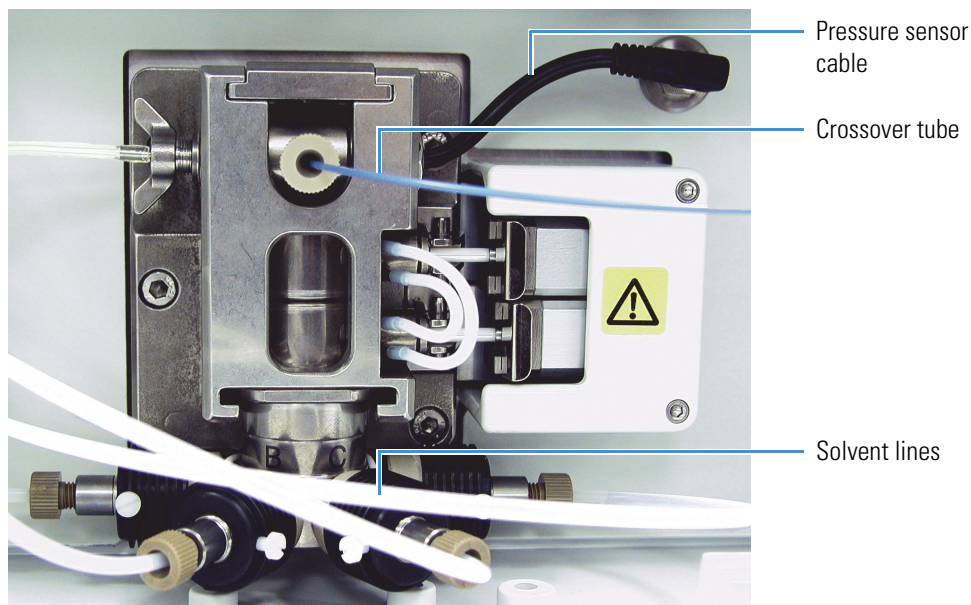
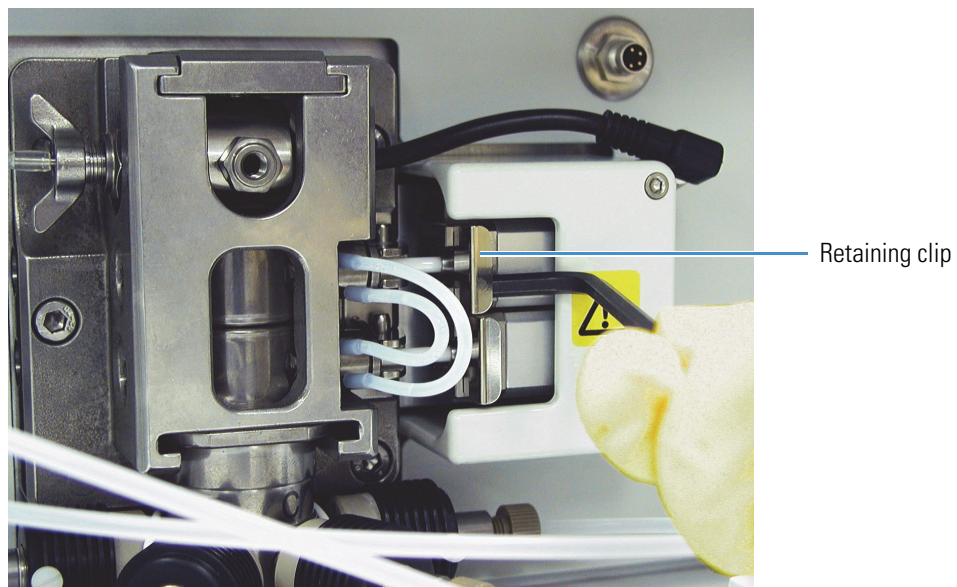


Figure 31. Using leverage to remove the piston spring clips



7. Detach the LDA from the pump by loosening the three captive screws securing the LDA body to the front of the pump. Carefully pull the LDA away from the pump by 8 to 10 cm (3 to 4 in.). See [Figure 32](#).

You cannot pull the LDA much farther away from the pump than this because the proportioning valve coils are connected to the pump by wires. See [Figure 33](#).



CAUTION Take care not to pinch the wires that connect the coils to the MS pump. Pinching a connecting wire can irreversibly damage the thin electrical wire housed within the white insulating coating. The damage might not be apparent until you attempt to run the MS pump.

Note Ideally your Surveyor MS Pump Plus is located on a clean, accessible benchtop where you can perform maintenance. If, however, your MS pump is located in an area in which it is inappropriate or difficult to perform maintenance, remove the LDA from the pump and place it on a clean, accessible benchtop before you disassemble it.

8. If necessary, remove the LDA from the pump. Use a flat-blade screwdriver to loosen the proportioning valve coils, and then pull the coils off the valves.

Figure 32. Location of the three captive screws

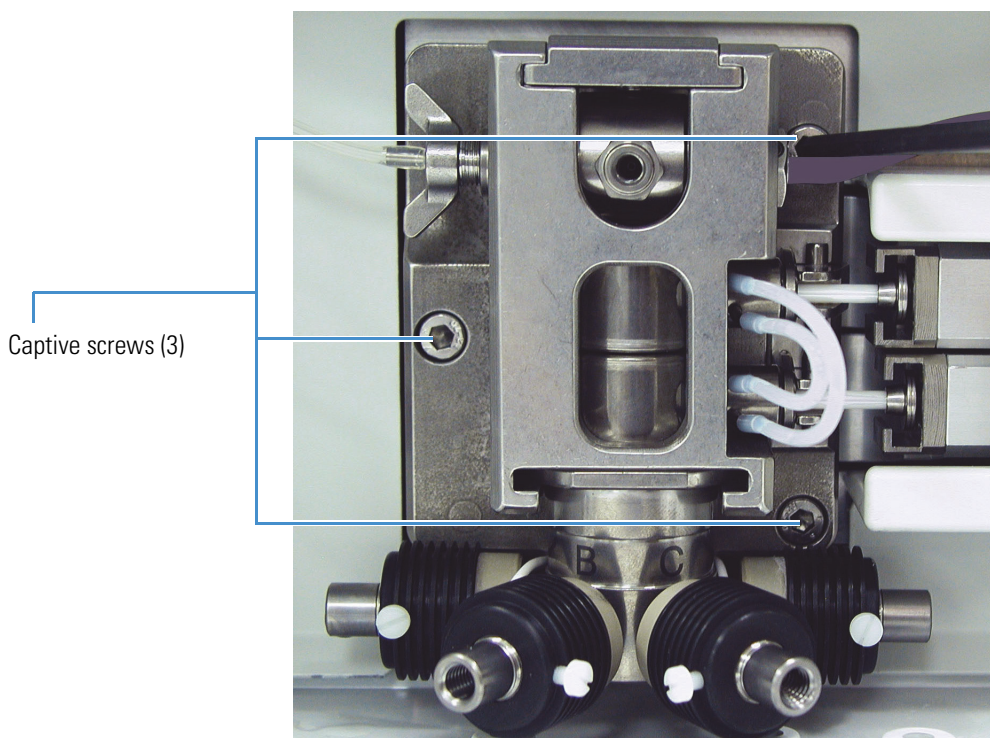
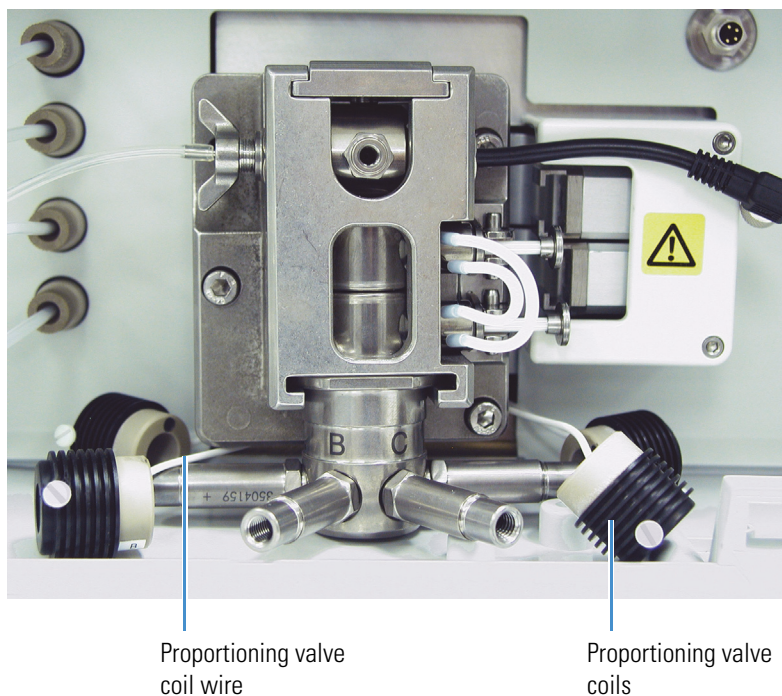


Figure 33. Removing the proportioning valve coils



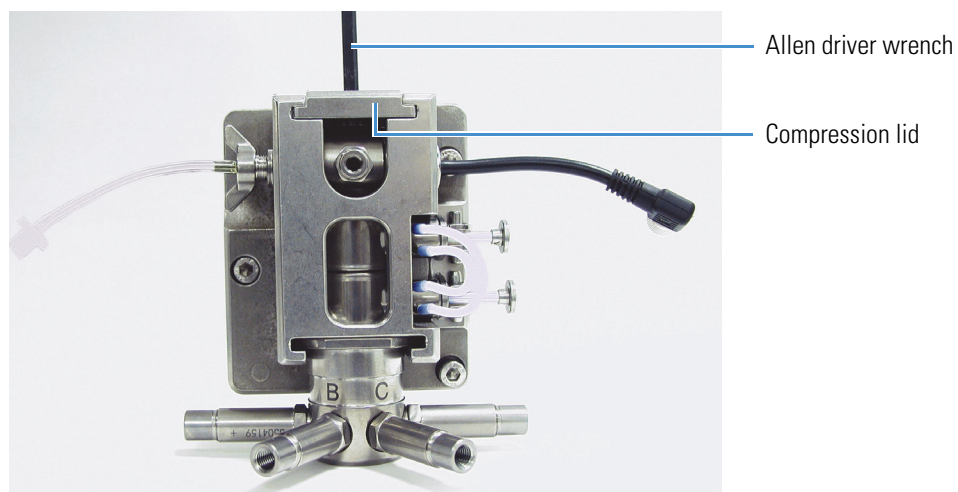
Disassembling the LDA

To prevent contamination of the internal surfaces of the LDA, make sure that you have a clean benchtop to work on before you disassemble the LDA.

❖ To disassemble the LDA

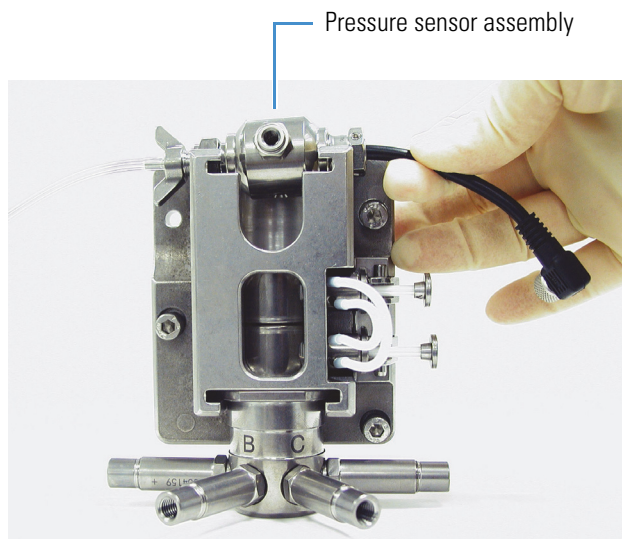
1. Use the 4 mm balldriver to loosen the top setscrew, and then slide the LDA compression lid out. See [Figure 34](#).

Figure 34. Removing the LDA compression lid



2. Remove the pressure sensor assembly by lifting it up and away from the LDA. See [Figure 35](#).

Figure 35. Removing the pressure sensor assembly



3. Remove the two pistons by pulling them to the right. See [Figure 36](#).
4. Use the 4 mm balldriver to loosen the fastening screws. The fastening screws are on the left side of the LDA and are attached to the holding hooks. See [Figure 37](#).

Figure 36. Removing the pistons from the LDA

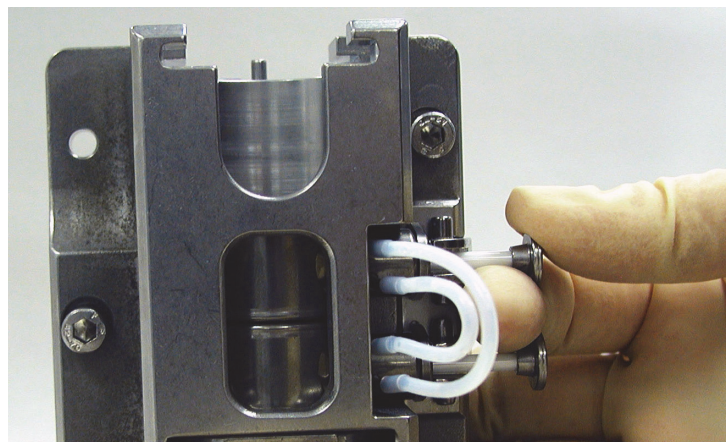
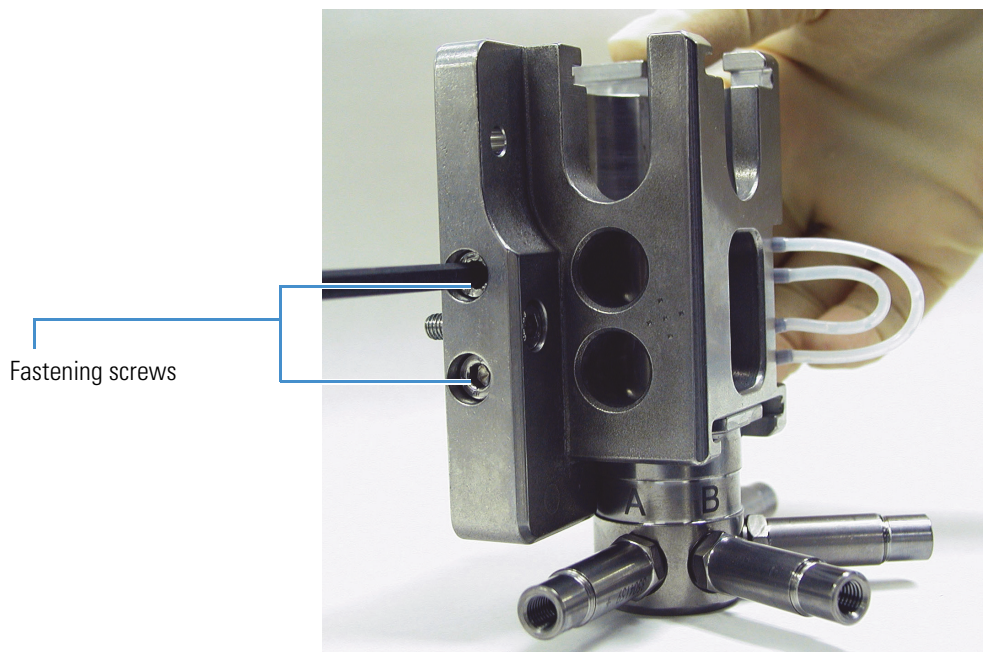
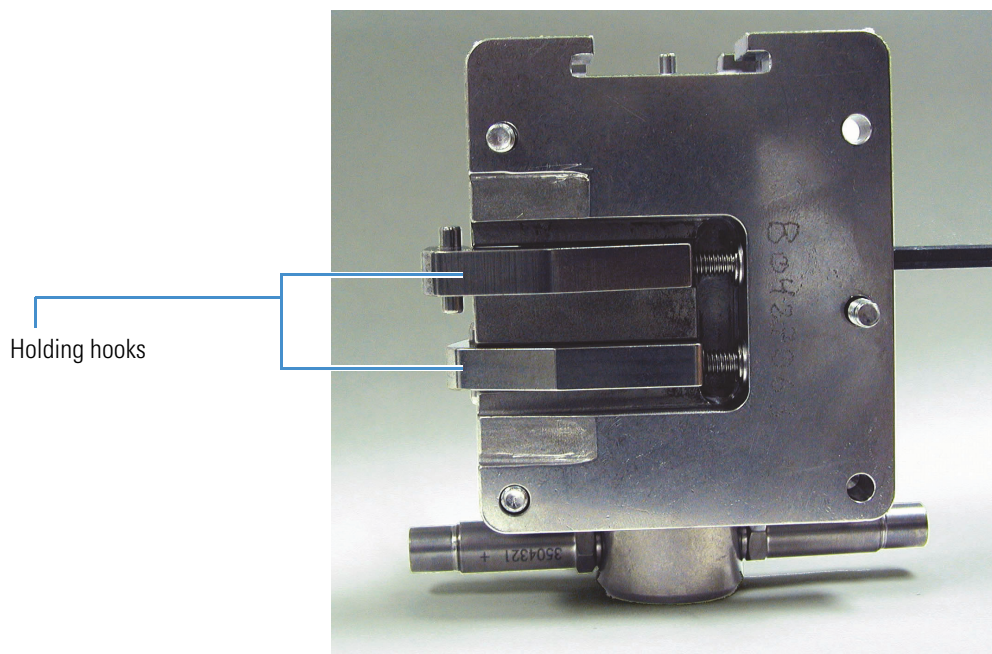


Figure 37. Loosening the two fastening screws



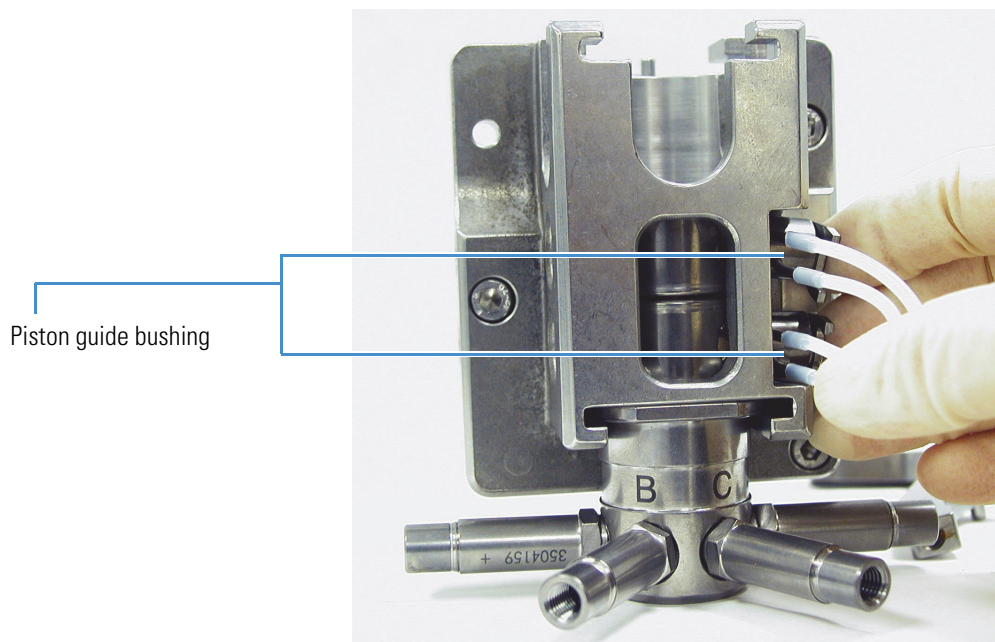
To provide enough clearance to remove the piston guide bushings, loosen the screws by at least six turns. As you loosen the screws, the holding hooks move to the right. If you prefer, you can detach the holding hooks completely with approximately nine turns on the screw fasteners. See [Figure 38](#).

Figure 38. View of loosened holding hooks



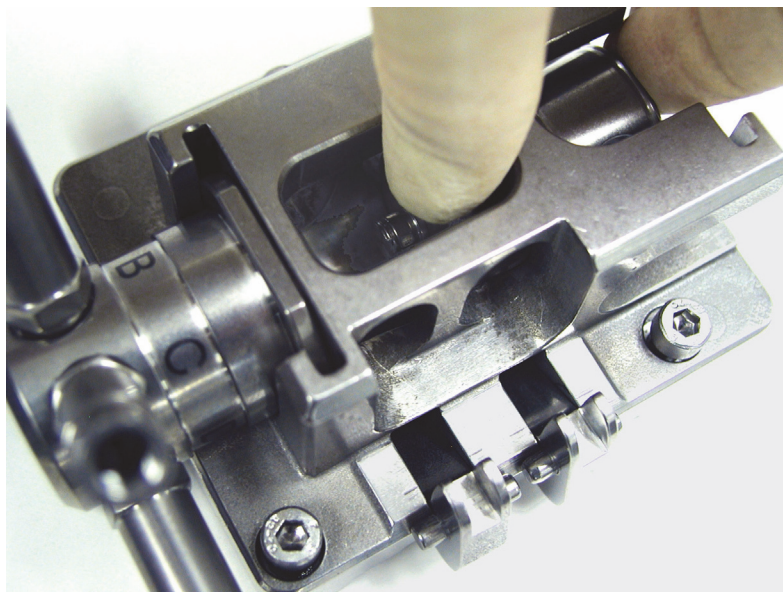
5. Remove the piston guide bushings and backing rings and place them on a clean surface. See [Figure 39](#).

Figure 39. Removing the piston guide bushings



6. Slide the two piston chambers out of the LDA body. To prevent canting, (tilting out of alignment, which can cause the piston chambers to get wedged within the inner cylinder of the LDA body) keep the piston chambers together as you remove them from the LDA body. See [Figure 40](#).

Figure 40. Removing the piston chambers



Cleaning the LDA Components

You can clean most LDA components without disassembling the LDA simply by running the pump for one hour with an appropriate solvent on all four channels (see “Cleaning the LDA Components without Disassembling the LDA” on page 38).

After you disassemble the LDA, concentrate on cleaning the following components:

- Check valves (see “Cleaning the Check Valves,” below)
- Pistons (see “Cleaning the Pistons” on page 48)
- Piston chambers and inlet module (see “Cleaning the Piston Chambers and the Inlet Module” on page 48)

These are the most likely components to have problems with contamination. The primary piston seals cannot be cleaned. If the primary piston seals contain particulate matter, replace them.

Cleaning the Check Valves

To clean the check valves, Thermo Fisher Scientific recommends using a 10 mL syringe with a very short piece of silicon tubing at the Luer exit. See Figure 41. With this syringe, gently flush the check valves with the following solvents: 5 mL acetone, followed by 5 mL of distilled water with a few drops of detergent, and then 5 mL methanol. If you cannot achieve free flow through the check valve, replace it.



CAUTION Thermo Fisher Scientific does not recommend cleaning check valves with an ultrasonic bath, because this procedure does not always extract fibrous particles from the check valves. You might even damage the small ruby balls inside by using an ultrasonic bath.

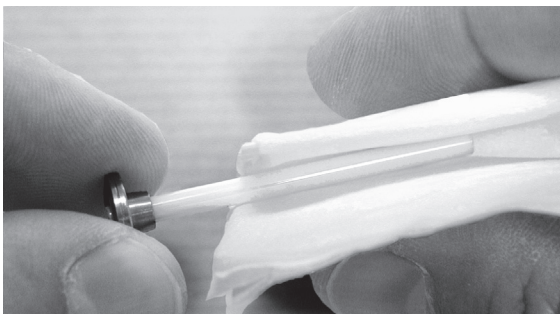
Figure 41. Cleaning the check valve



Cleaning the Pistons

If precipitate is visible on the pistons (P/N 00201-11324), clean them with a soft tissue and cleaning paste. See [Figure 42](#).

Figure 42. Cleaning a piston



Cleaning the Piston Chambers and the Inlet Module

Clean the piston chambers if the primary piston seals have started to decompose and you notice black, fibrous, particles in the chamber.

❖ To clean the piston chambers and the inlet module

1. Insert the seal removal tool into the used primary piston seal. Use a twisting motion to pry the seal out of the piston chamber. See [Figure 43](#). The primary piston seals cannot be reused. After you remove them from the piston chamber, throw them away.

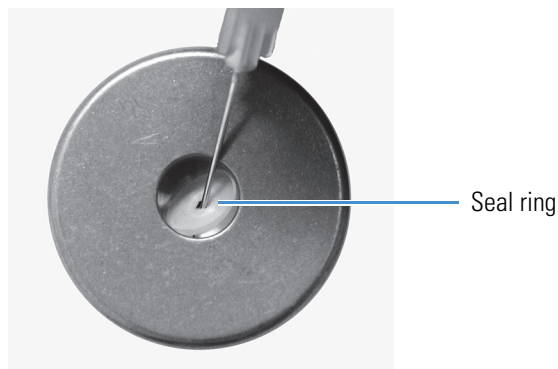
Figure 43. Removing the used primary piston seal



Note Each side of the piston chambers and the top bore of the inlet module contain a check valve seal ring for a total of five seal rings. Inspect the seal rings inside the top and the bottom of the piston chambers.

2. Use a needle to remove the check valve seal rings (P/N 00107-18118) as shown in [Figure 44](#). Be careful not to scratch the piston chamber with the needle as you remove its seal ring, as scratching the piston chamber might cause leaking.

Figure 44. Using a needle to remove the check valve seal ring



3. Place the piston chambers and the inlet module in an ultrasonic bath. First use water (to remove any buffer precipitate) and then use acetone.
4. After cleaning the piston chambers and the inlet module, reinstall the seal rings if they do not show signs of damage. Unlike the seal rings, the primary piston seals cannot be reused. Replace the primary piston seals before reassembling the LDA.

Go on to the next topic: [Replacing the Primary Piston Seals](#).

Replacing the Primary Piston Seals

You must replace the primary piston seals when you disassemble the LDA.

❖ To replace the primary seals

1. Soak a pair of new primary piston seals (springs up) for five minutes in methanol. Soaking the seals removes air from the springs.

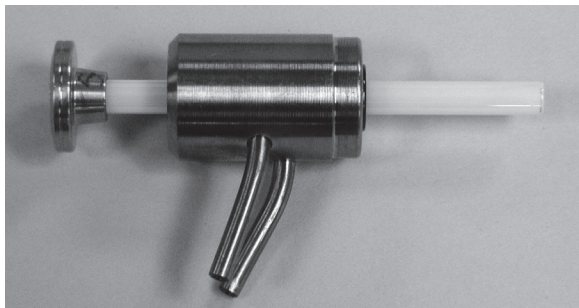


CAUTION Do not place the seals in an ultrasonic bath. Placing the seals in an ultrasonic bath can ruin the smooth surface of the seals.

2. If you have not already done so, do the following:
 - a. Remove the LDA from the pump (see [“Removing the LDA from the Pump”](#) on [page 40](#)).
 - b. Disassemble the LDA (see [“Disassembling the LDA”](#) on [page 43](#)).
 - c. Remove the used piston seal (see [“Cleaning the Piston Chambers and the Inlet Module”](#) on [page 48](#)).

3. Install a new primary piston seal (P/N 00107-18110):
 - a. Create a seal insertion tool by inserting a clean piston into a piston guide bushing as shown in [Figure 45](#).

Figure 45. Seal installation tool (piston and piston guide bushing)



- b. Slide the new seal onto the piston. Be sure that the spring is facing out, so that it faces the piston chamber when the seal is installed in the chamber. See [Figure 46](#).

Figure 46. View of the spring face



- c. Insert the piston, which is holding the seal, into the piston bore.
 - d. Push the end of the piston guide bushing into the piston bore to make sure that the seal is properly seated in the piston chamber. See [Figure 47](#) and [Figure 48](#).
4. Repeat steps 2 and 3 with the second piston chamber.

Figure 47. Pushing the piston guide bushing into the piston bore

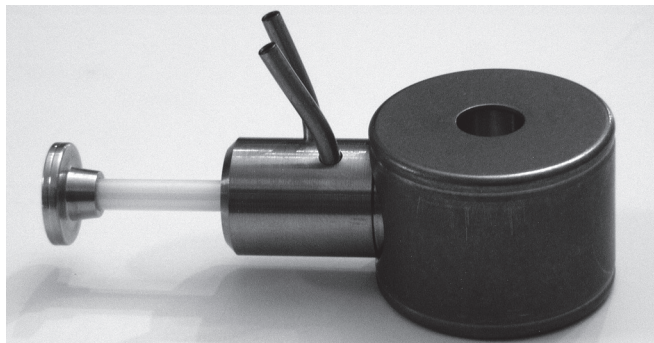
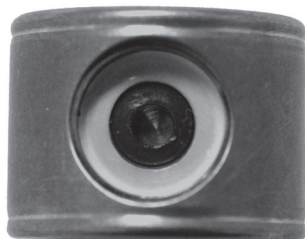


Figure 48. View showing the piston bore with a properly seated seal



Reassembling the LDA

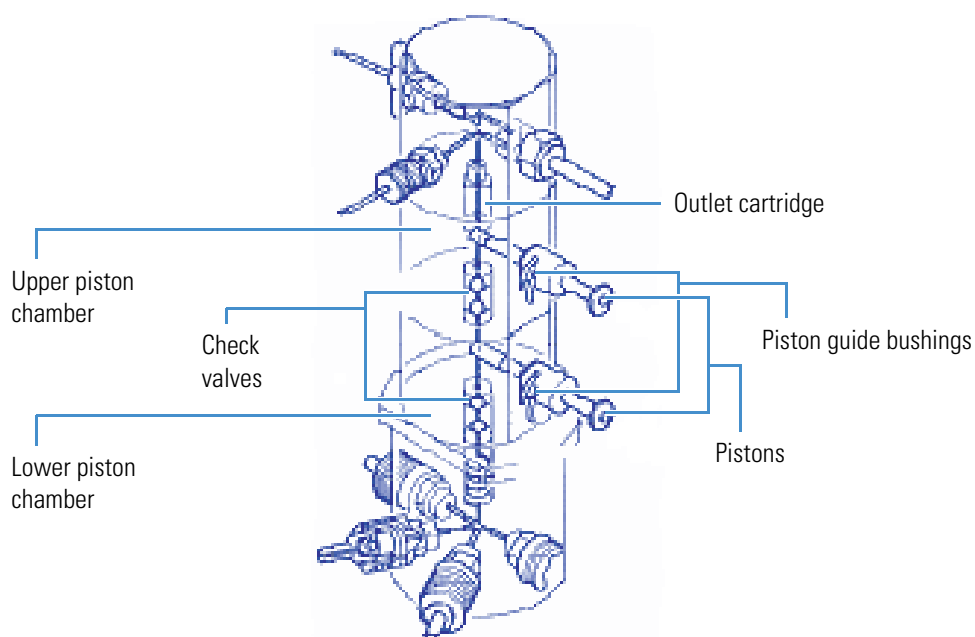
The LDA is reassembled in three stages. First, the stacked components that fit within the LDA body are reinstalled. Next, the piston guide bushings, the backing rings, and the pistons are reinstalled. And finally, the compression lid and the left side fasteners are tightened with the 4 mm Allen wrench to prevent leaking.

❖ **To reassemble the LDA, follow these steps**

- [Step 1: Reinstalling the Stacked Components of the LDA](#)
- [Step 2: Reinstalling the Piston Guide Bushings and the Pistons](#)
- [Step 3: Tightening the LDA to Prevent Leakage](#)

If they are mounted correctly, the parts of the LDA form a perfect stack to ensure a smooth piston movement. A schematic diagram of the LDA is shown in [Figure 49](#).

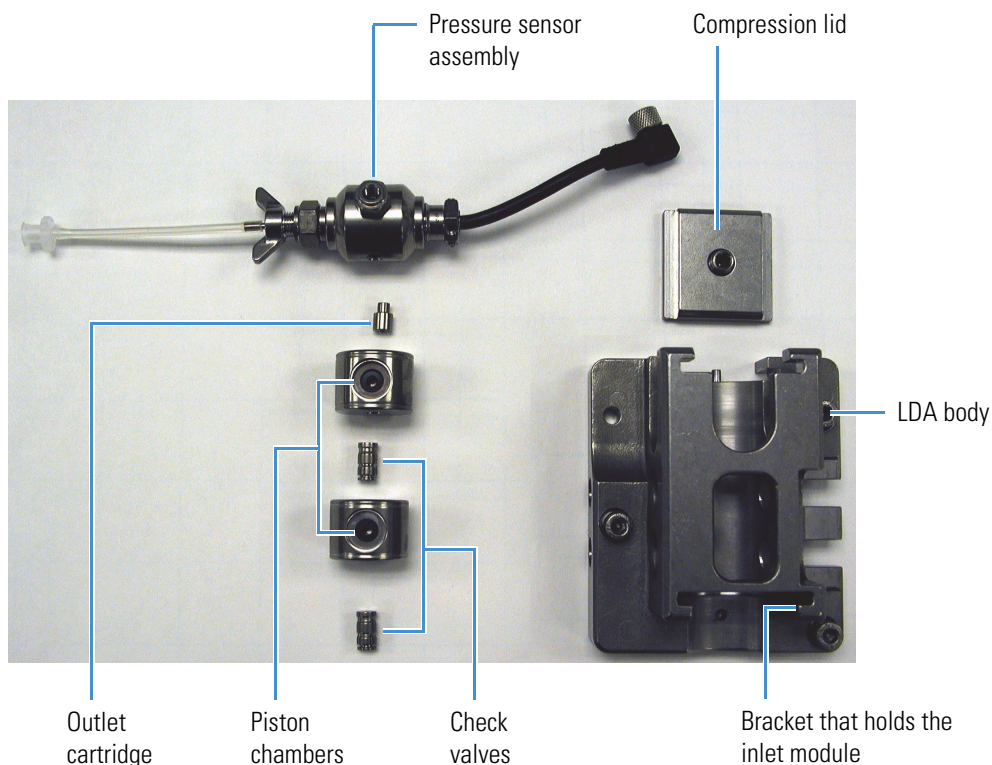
Figure 49. LDA schematic diagram



Step 1: Reinstalling the Stacked Components of the LDA

The stacked components of the LDA form a linear flow path from the inlet module on the bottom of the stack to the pressure sensor assembly on the top of the stack. See [Figure 50](#).

Figure 50. LDA stack components



❖ **To reinstall the components of the stack**

1. Take a check valve (P/N 00110-05110), noting the flow direction (flow is from bottom to top of the LDA), and drop it into place in the inlet module. See [Figure 51](#) and [Figure 52](#).

Figure 51. Check valve, showing the flow direction

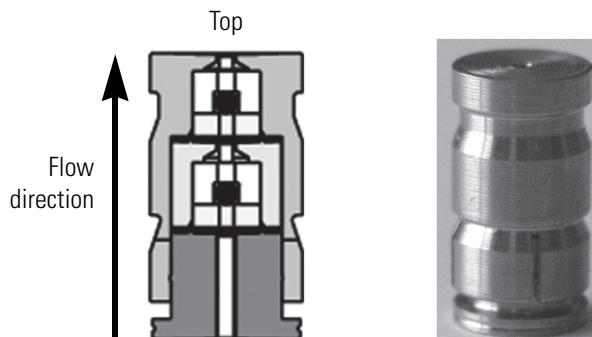
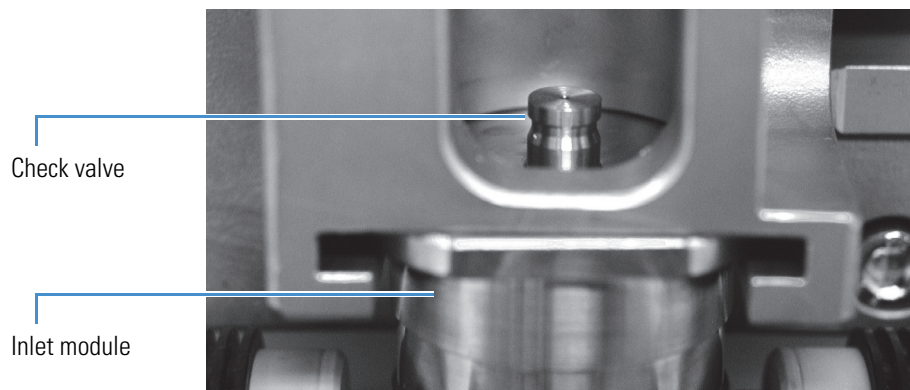
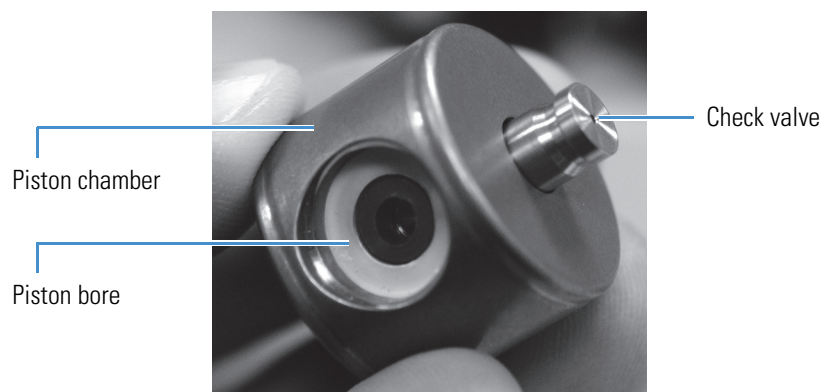


Figure 52. Inserting the check valve into the inlet module



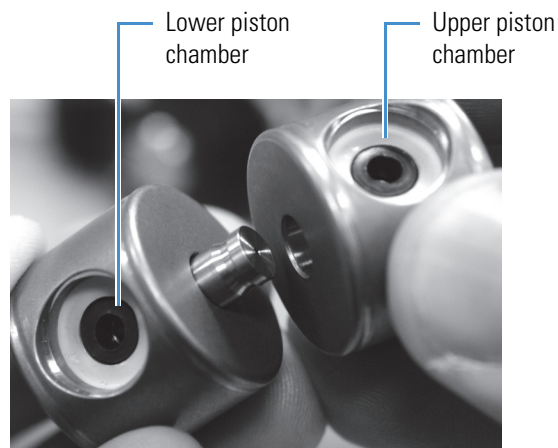
2. To prevent canting of the chambers as you insert them into the LDA body, create a stack of the two piston chambers with the remaining check valve in the middle:
 - a. Drop the second check valve into the first piston chamber (noting flow goes up as [Figure 51](#) shows). See [Figure 53](#).

Figure 53. Inserting check valve into piston chamber



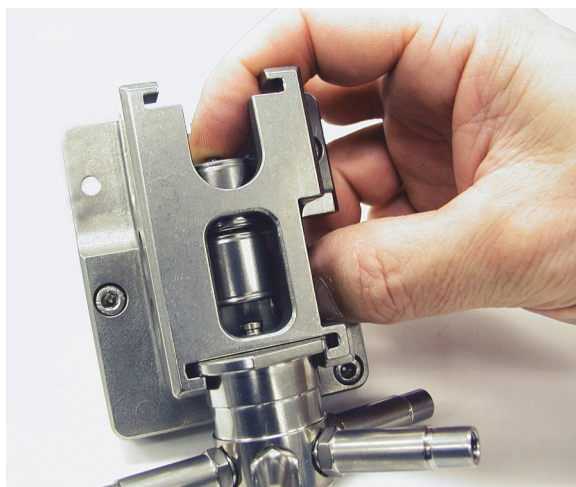
- b. Take the second chamber and slide it on top. Orient the second chamber so that its piston bore is aligned with the piston bore on the first chamber. See [Figure 54](#).

Figure 54. Aligning the piston chambers



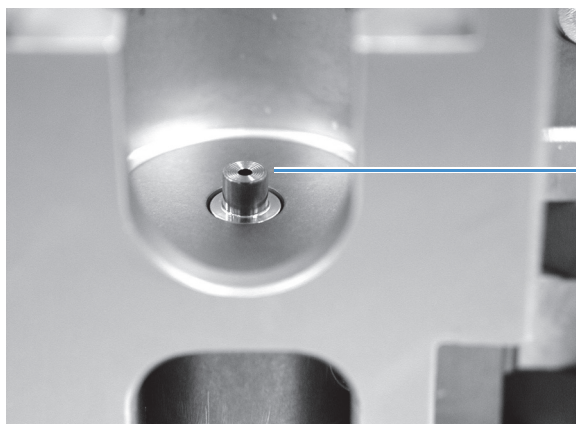
3. Slide the stacked piston chambers into the LDA body so that the piston bores are to the right. See [Figure 55](#). Ensure that the stack is oriented with the direction of flow from bottom to top.

Figure 55. Installing the piston chambers



4. Insert the outlet cartridge, with the narrow end up, into the upper piston chamber. See [Figure 56](#).

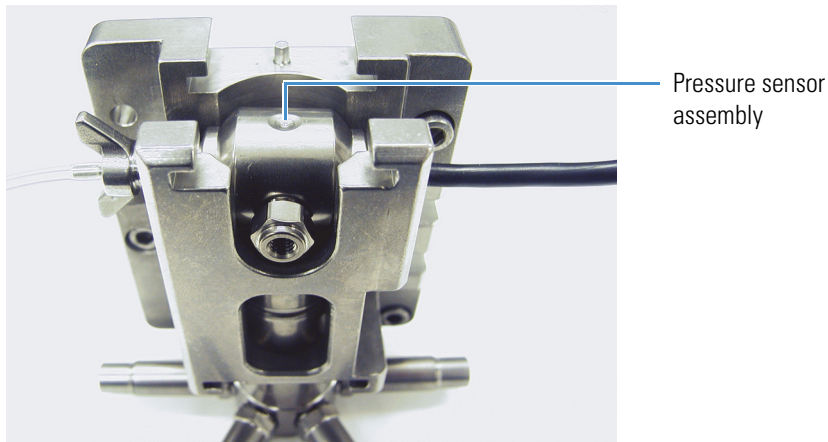
Figure 56. Reinstalling the outlet cartridge



Outlet cartridge

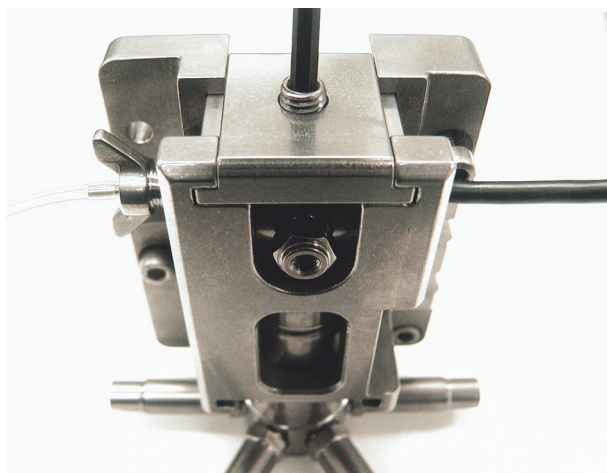
5. Align the bore on the bottom of the pressure sensor assembly with the outlet cartridge, and then set the pressure sensor assembly in place. See [Figure 57](#).

Figure 57. Installing the pressure sensor



6. Slide the compression lid back into place. Tighten the screw until it is snug and then loosen it by half a turn. You tighten the screw later once the pistons and the piston guide bushings are in place. See [Figure 58](#).

Figure 58. Tightening the compression lid



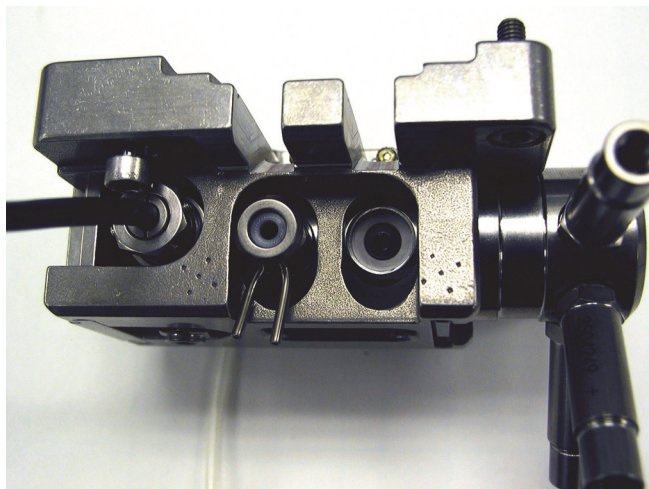
Step 2: Reinstalling the Piston Guide Bushings and the Pistons

After you have reinstalled the stacked components of the LDA, you are ready to reinstall the piston guide bushings and the pistons.

❖ To install the guide bushings and the pistons

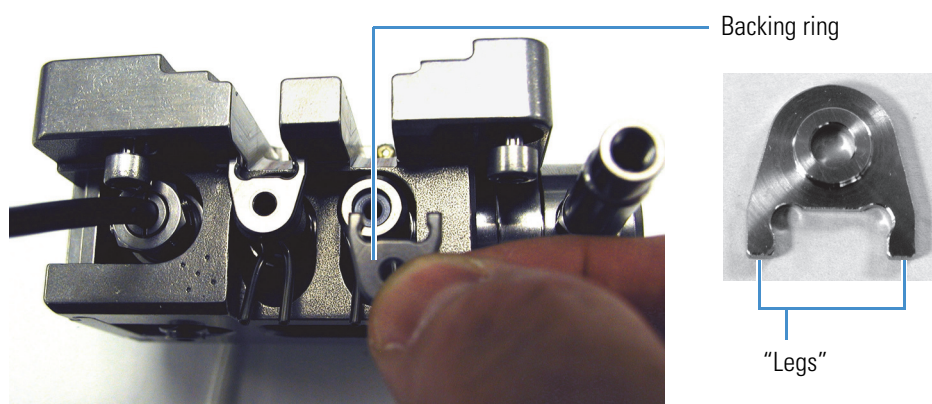
1. Insert the piston guide bushings into the piston bores. The end of the piston guide bushing that contains the red ruby bearing fits into the piston bore. See [Figure 59](#).

Figure 59. Inserting the piston guide bushing into the piston bore



2. Insert the two A-shaped backing rings into the other end of the piston guide bushings. Make sure that the “legs” of the backing rings face toward the holding hooks as shown in [Figure 60](#).

Figure 60. Installing the backing rings

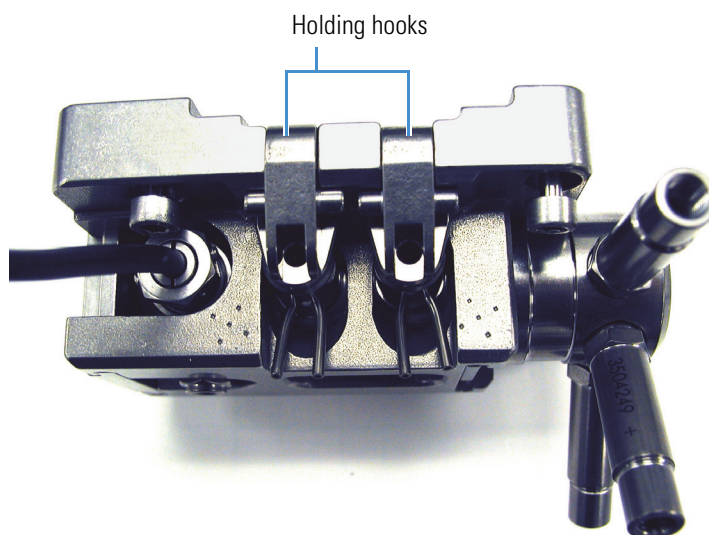


3. Alternately tighten the fasteners on the left side of the LDA until the holding hooks are close enough to prevent slippage of the piston guide bushings and the backing rings. See [Figure 61](#) and [Figure 62](#).

Figure 61. Tightening the left-side fasteners, as viewed from the back of the LDA body

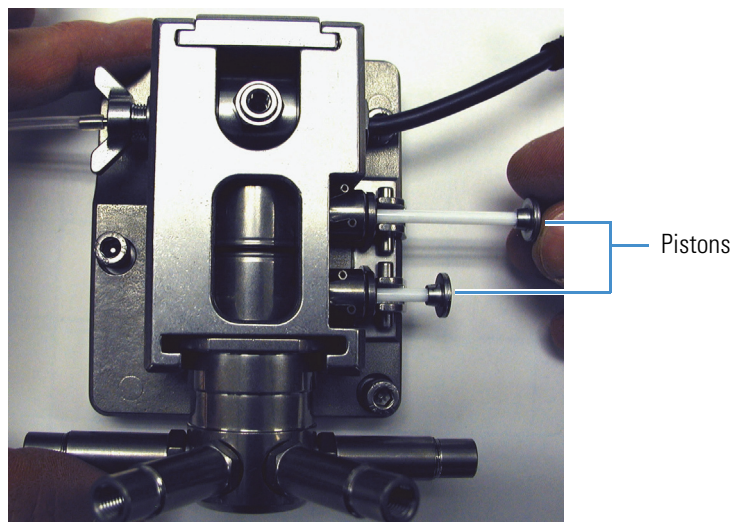


Figure 62. View of piston guide bushings and backing rings held snugly by holding hooks



4. Insert the two pistons into the holes in the backing rings. Continue to push them into the piston bores until they contact the bottom. As you push the pistons into the bores, you can feel interference due to the seals. If the pistons appear uneven, loosen the left side fasteners and readjust the pistons until they contact the bottom of the piston bores. See [Figure 63](#).

Figure 63. Installing the pistons



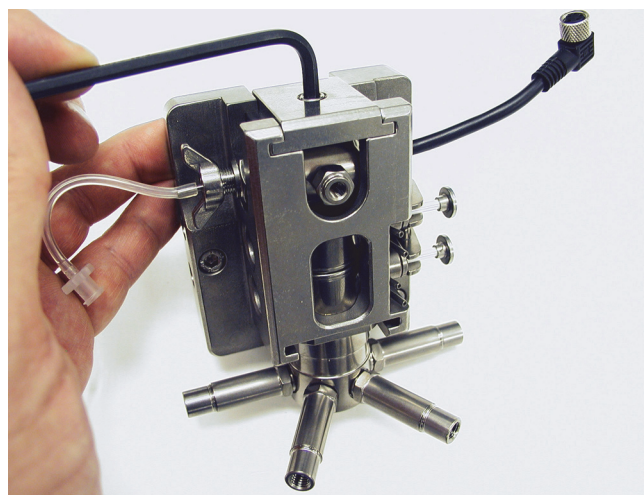
Step 3: Tightening the LDA to Prevent Leakage

After you have reinstalled all the components of the LDA, you are ready to perform the final tightening that prevents leakage.

❖ **To tighten the compression lid and the left side fasteners**

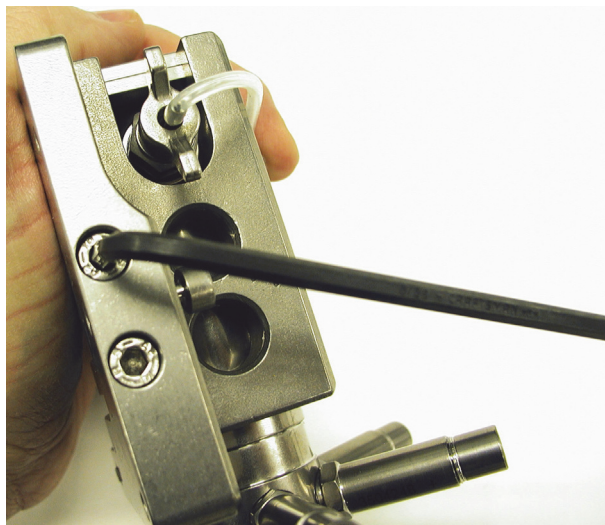
1. Using an Allen wrench, tighten the compression lid by an additional quarter to half turn, and then retest the piston movement. See [Figure 64](#).

Figure 64. Final tightening of the compression lid



2. Using an Allen wrench, tighten the two left side fasteners by an additional quarter to half turn, and then verify that the pistons move in and out easily, but with some interference due to the seals. See [Figure 65](#).

Figure 65. Final tightening of the left-side fasteners



Reinstalling the LDA

After you reassemble the LDA, you are ready to reinstall the LDA onto the MS pump.

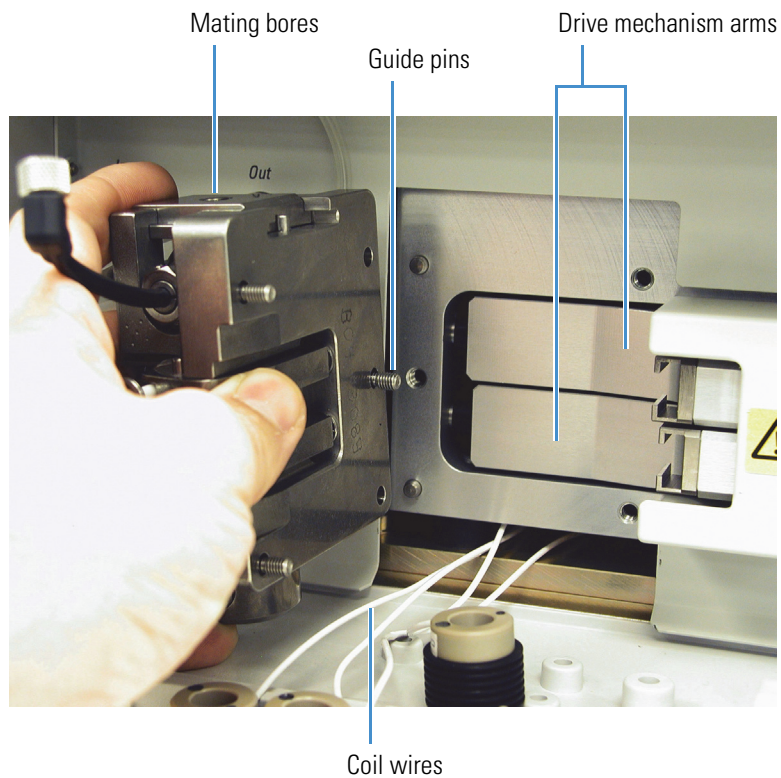
❖ To reinstall the LDA

1. Reattach the coils to the proportioning valves.
2. Position the LDA onto the front of the pump. The two guide pins on the pump drive align with the two mating bores on the LDA housing. Before you attach the LDA to the pump, make sure that the wires connected to the proportioning valve coils are located in the slot below the arm drive mechanism. See [Figure 66](#).



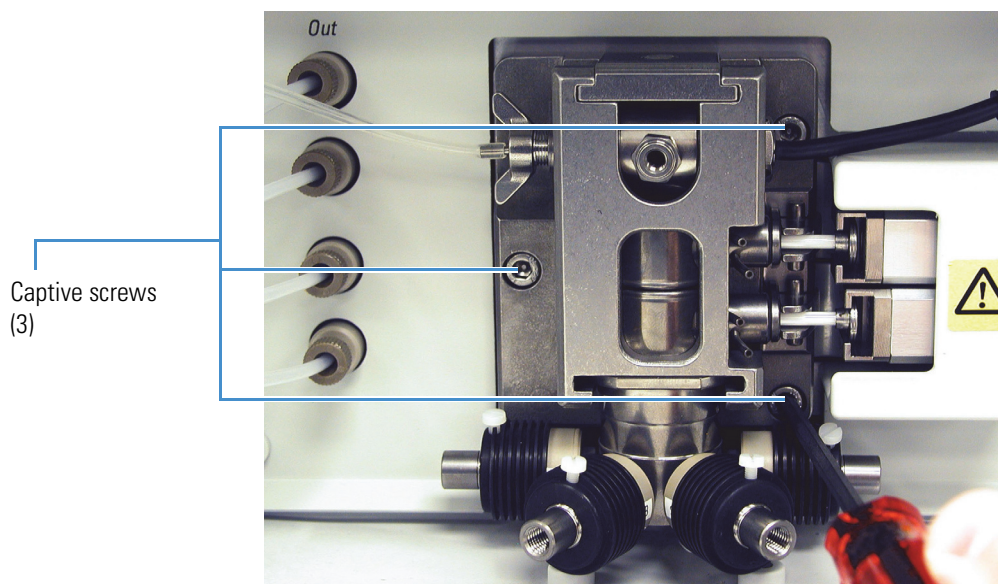
CAUTION Although the damage might not be visible, pinching one of these connecting wires can irreversibly damage the internal copper wire, which conducts electricity to magnetize the coil. This damage causes the proportioning valve to malfunction.

Figure 66. Positioning the LDA on the pump



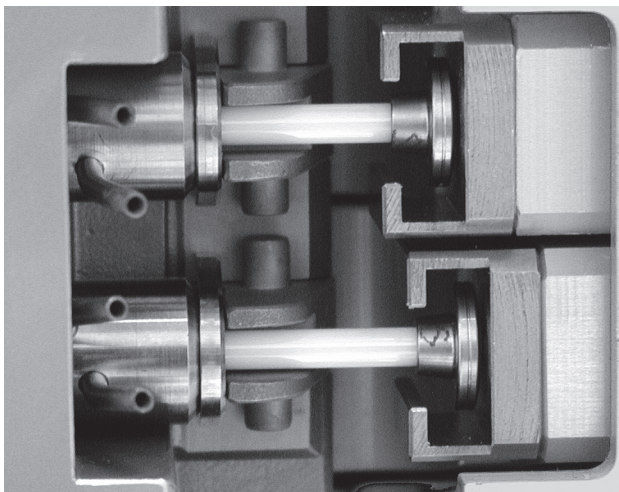
3. Alternately tighten the three captive screws that hold the LDA body to the pump. See [Figure 67](#).

Figure 67. Attaching the LDA to the pump



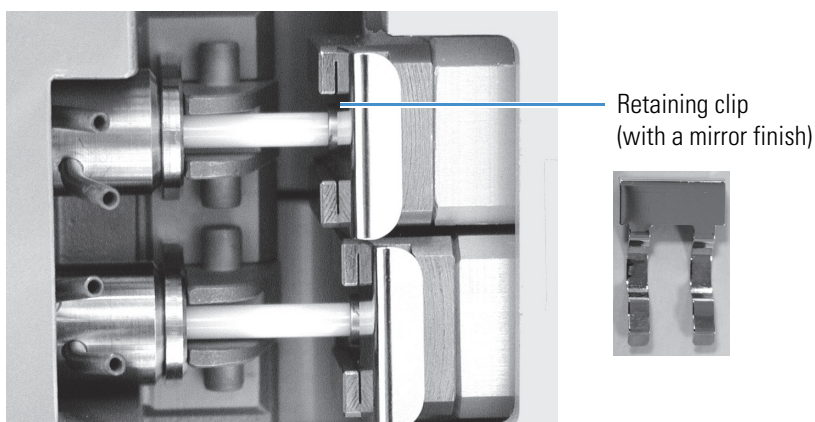
4. Push the piston heads toward the arms of the drive mechanism. Leave a gap for the piston retaining clips between the piston head and the brackets at the ends of the arms. See [Figure 68](#).

Figure 68. Pushing the piston heads toward the drive mechanism arms



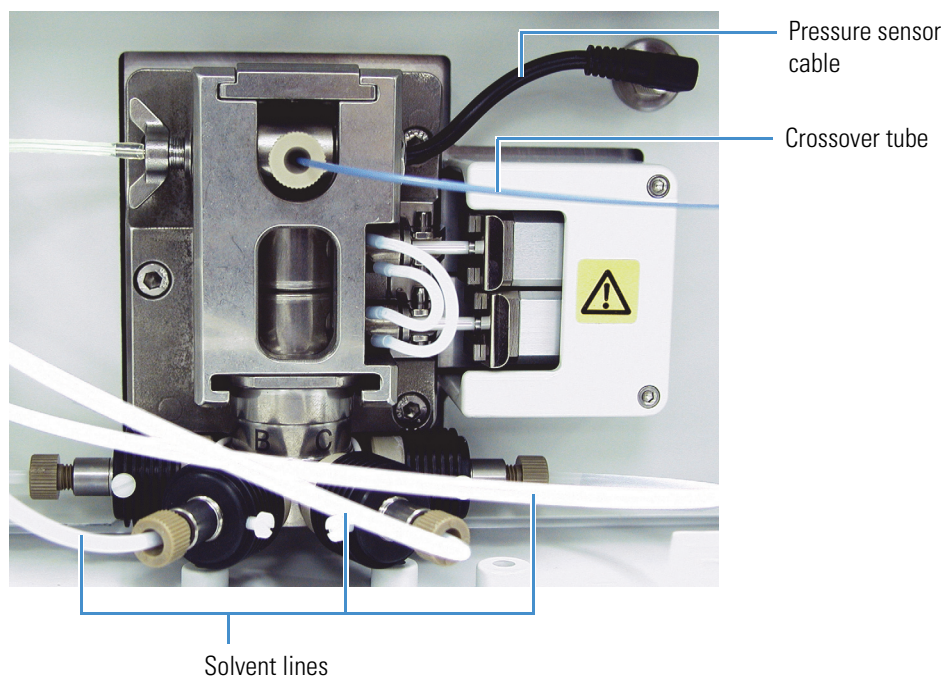
5. Insert the two piston retaining clips between the piston heads and the brackets of the drive mechanism arms. See [Figure 69](#).

Figure 69. Installing the piston retaining clips



6. Reconnect the following:
 - Connect the pressure sensor cable to the pump (see [Figure 70](#)).
 - Connect the crossover tube to the pressure sensor assembly (see [Figure 70](#)).
 - Connect the solvent lines to the four proportioning valves (see [Figure 70](#)).
7. Turn on the solvent flow and check for leaks.
8. Perform the pump pulsation and proportioning tests described in [Chapter 3, "Diagnostics,"](#) to ensure that your pump is functioning properly.

Figure 70. Completed assembly

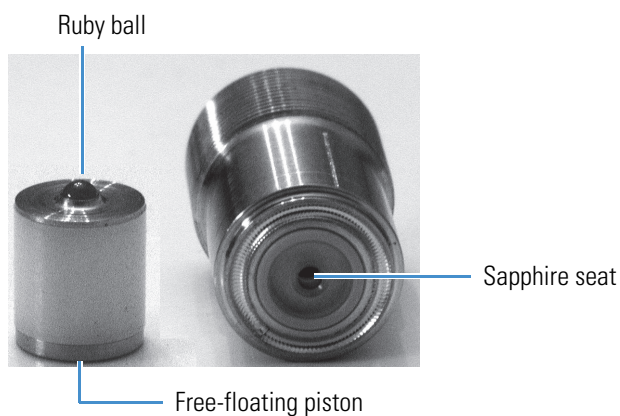


Servicing the Proportioning Valves

Each proportioning valve of the Surveyor MS Pump Plus contains an electromechanically actuated ball and seat valve. The ruby ball portion of the valve sits on one end of a magnetic piston that floats within the body of the valve. The sapphire seat, against which the ball rests, is located on the backside of the valve face. See [Figure 71](#).

The ruby ball normally rests against the sapphire seat, shutting off the solvent flow. Electrical pulses from the coils surrounding the valve cause the magnetic piston to glide back and forth within the body of the valve, allowing solvent to flow through the valve and into the inlet module.

Figure 71. Ruby ball and sapphire seat



Occasionally, after extended periods of disuse, the free-floating piston inside the proportioning valve freezes against its sapphire seat. When this occurs, dislodging the frozen valve requires a greater mechanical force than the magnetic field applied by the actuating coils.

You can dislodge the free-floating piston inside the valve with the proportioning valve tool supplied in the accessory kit. See [Figure 72](#). This tool fits through the solvent inlet in the end of the valve face. Push the tool (which is a piece of stainless steel wire) into the valve until you contact the piston, and then gently push the piston away from the valve face.

Figure 72. Proportioning valve, shown with the valve tool



Replacing the Solvent Reservoir Filter

The solvent reservoir filter removes particulate matter from the mobile phase before it enters the solvent line. Eventually particulate matter from the mobile phase builds up on the filter and restricts the flow of solvent. A clogged solvent reservoir filter can cause the formation of air bubbles in the solvent line, erratic flow rates, and an abnormally low system backpressure.

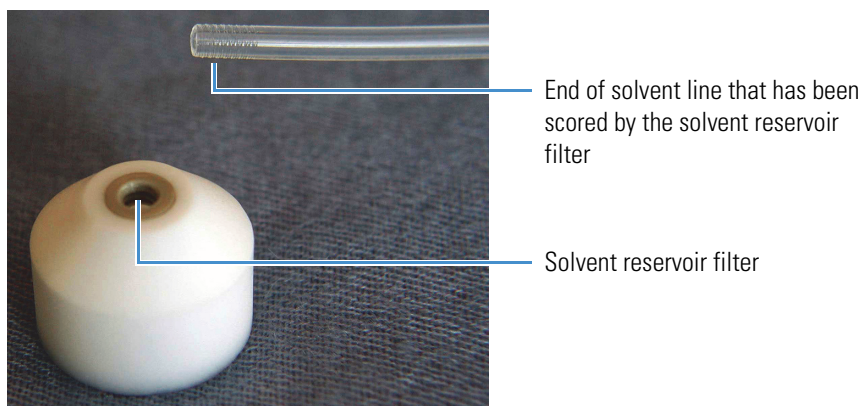
❖ To determine if the solvent reservoir filter is clogged

1. Remove the filter from the solvent line
2. Monitor the system backpressure as the pump draws solvent into the solvent line.

If the pressure trace is normal in the absence of the solvent reservoir filter, replace the filter.

Thermo Fisher Scientific offers two solvent reservoir filters: a 10 micron pore size Teflon filter (P/N A4258-010) and a 10 micron pore size stainless steel filter (P/N A4929-010). [Figure 73](#) shows the Teflon® filter. The Teflon filter screws onto the end of the solvent line.

Figure 73. Solvent reservoir filter and solvent line



Troubleshooting

If properly maintained, the Surveyor MS Pump Plus can operate free of trouble for many years. Most pump problems can be avoided by simple, periodic maintenance, as described in [Chapter 4, “Maintenance.”](#) However, in the event that an error message is displayed or if a mechanical or electrical failure is suspected, you might be able to diagnose the source of the problem. If the diagnosis indicates that a problem exists with serviceable parts inside the pump, contact a Thermo Fisher Scientific service representative.

Contents

- [Troubleshooting Tips](#)
- [General System Troubleshooting](#)
- [Surveyor MS Pump Troubleshooting Guide](#)
- [Pump Pulsation](#)
- [Baseline Noise at Low Pressures \(under 40 Bar\)](#)
- [Insufficient Solvent Flow from the Pump Outlet](#)
- [Irreproducible Retention Times](#)

For more detailed chromatographic troubleshooting, see any HPLC troubleshooting reference book or call your Thermo Fisher Scientific service representative.

Troubleshooting Tips

Begin the diagnosis of the problem by performing the following steps:

Eliminate all other possible sources of trouble. Before you invest time trying to diagnose a suspected pump problem, you should verify that the pump is the only source of difficulty. Systematically eliminate all other instruments in your Surveyor HPLC system as the source of trouble. If you are not sure which component of your chromatography system is responsible for poor system performance, see [“General System Troubleshooting”](#) on [page 66](#) for some useful suggestions.

When you are sure that it is the pump. When you have isolated the pump as the only remaining source of difficulty, proceed to the troubleshooting table for help in determining the exact cause of the problem.

General System Troubleshooting

The following section can be helpful in the initial investigation of a problem with the Surveyor Plus HPLC system.

❖ **No flow**

1. Check mobile phase connections.
2. Check for leaks.
3. See problem 8 on [page 69](#).

❖ **High Backpressure**

1. Check flow rate and system/column specifications.
2. Check for tubing or column blockage.
3. Check for blockage in the injection valve.
4. See “[Surveyor MS Pump Troubleshooting Guide](#)” on [page 68](#).

❖ **Unstable Baseline or Drift**

1. System column not equilibrated; allow more equilibration time.
2. Refer to the troubleshooting guide for your detector.
3. See problem 7 on [page 69](#).

❖ **Baseline Noise**

1. Check for air bubbles in the system.
2. Check for system solvent contamination.
3. The pulse dampener is contaminated; prime the pulse dampener.

❖ **No Peaks**

1. Check the detector and data system connections.
2. Refer to the troubleshooting guide for your autosampler.
3. Check sample retention with chromatographic conditions.
4. See problem 8 on [page 69](#).

❖ **Contaminating / Ghost Peaks**

1. Clean the system and the column.
2. Refer to the troubleshooting guide for your autosampler.

❖ **Poor Peak Shape**

1. Check the system for leaks.
2. Check the fittings and tubing lengths.
3. Check the column performance.
4. Refer to the troubleshooting guide for your autosampler.
5. Refer to the troubleshooting guide for your detector.

❖ **Poor Retention Time Reproducibility**

1. Check the system for leaks and bubbles.
2. System/column not equilibrated and more time should be allowed.
3. Check the column performance.
4. Refer to the troubleshooting guide for your detector.
5. See problem 7 on page [page 69](#).

❖ **Poor Peak Area Reproducibility**

1. Check the column performance.
2. Refer to the troubleshooting guide for your autosampler.

❖ **No Instrument or Device Control**

1. Check the cable connections.
2. Check the system configuration.
3. Refer to the troubleshooting guide for each instrument.

Surveyor MS Pump Troubleshooting Guide

The information in [Table 4](#) can help you to identify specific pump-related problems.

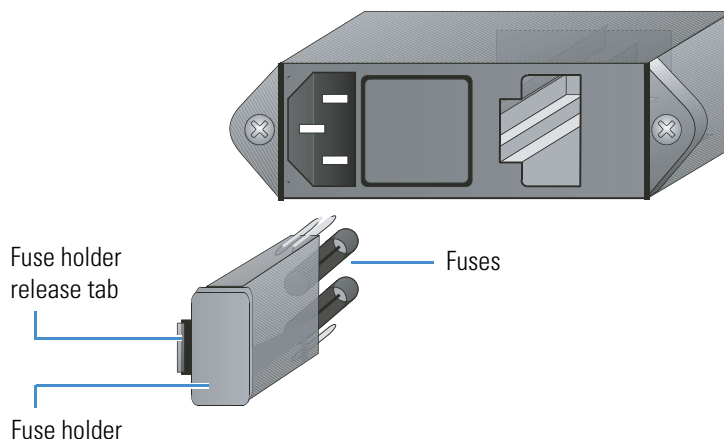
Table 4. Troubleshooting pump related problems (Sheet 1 of 2)

Problem	Cause	Diagnostic/Remedy
1. No response when power is switched on	a. Power cord not firmly installed	a. Reseat power cord
	b. Power cord defective	b. Replace cord
	c. Power entry module fuse is blown	c. Replace fuse (See Figure 74)
2. LEDs not illuminated	a. LED PCB defective	a. Contact Thermo Fisher Scientific service representative
	b. Loose LED cable	b. Reseat LED connector
3. Pump motor does not run	a. Max pressure set to zero	a. Change max pressure in Instrument Setup and/or Direct Control
	b. Motor defective	b. Contact Thermo Fisher Scientific service representative
	c. Motor cable disconnected	c. Contact Thermo Fisher Scientific service representative
	d. Defective drive circuit	d. Contact Thermo Fisher Scientific service representative
	e. Internal fuse blown	e. Contact Thermo Fisher Scientific service representative
4. Incorrect solvent composition	a. Proportioning valve dirty	a. Clean with acetic acid
	b. Proportioning valve cable loose	b. Reattach proportioning valve cable
	c. Proportioning valve timing not properly set	c. Contact Thermo Fisher Scientific service representative
	d. Proportioning valve defective	d. Contact Thermo Fisher Scientific service representative
	e. Proportioning valve drive circuit defective	e. Contact Thermo Fisher Scientific service representative
5. Pump does not start or stop remotely	a. Incorrect connection of synchronization cable	a. Reconnect synchronization cable

Table 4. Troubleshooting pump related problems (Sheet 2 of 2)

Problem	Cause	Diagnostic/Remedy
6. Proportioning valves do not click	a. Proportioning valve cable loose	a. Reattach proportioning valve cable
	b. Proportioning valve defective	b. Contact Thermo Fisher Scientific service representative
	c. Proportioning valve drive circuit defective	c. Contact Thermo Fisher Scientific service representative
	d. Proportioning valves are stuck	d. See “Servicing the Proportioning Valves” on page 62
7. Erratic flow rate	a. Restricted inlet tubing	a. Replace inlet tubing
	b. Restricted outlet tubing	b. Replace outlet tubing
	c. Dirty proportioning valve	c. Clean with acetic acid or replace proportioning valves
	d. Dirty check valve	d. Clean with acetic acid or replace check valves
	e. Leaking piston seal	e. Replace piston seals
8. No flow	a. Loose or disconnected solvent lines	a. Check for leaks in the solvent lines at the degassing inlets and outlets, the proportioning valve inlets, and the crossover tube connections
	b. Excessive air in the solvent lines	b. Remove the air from the solvent lines
	c. Open venting valve	c. Close the venting valve on the left side of the LDA

Figure 74. Power Entry Module



Pump Pulsation

Pump pulsation is the variation in pressure with time due to the motion of the pump pistons.

This section contains the following topics:

- [Normal Pressure Pulsation](#)
- [Abnormal Pressure Pulsation](#)

Normal Pressure Pulsation

The pump continuously monitors the system pressure and bases its pressure readout on a 30 s interval. Under normal operating conditions, the Surveyor MS Pump Plus shows very small variation in pressure. At pressures greater than 100 bar (1450 psi), the pump pulsation should be less than 1%.

The pressure status readback for the MS pump in the Information View is in standard deviation percent. Therefore, at backpressures greater than 100 bar, a value of 1.0 or less is acceptable. See “[Monitoring Pump Pulsation](#)” on [page 29](#) for instructions on monitoring the backpressure of your system from the data system.

Note Because the Surveyor MS Pump Plus does not use electronic damping, its pressure fluctuation often appears to be higher than that for conventional pumps.

Abnormal Pressure Pulsation

Abnormal pressure pulsations can normally be divided into two categories. Large pressure pulsations are generally considered to be those with a standard deviation in pressure of greater than 10 bar (147 psi). Minor pressure pulsations are those with a standard deviation in pressure of less than 10 bar.

Pressure Pulsations with SD over 10 Bar

Pressure pulsations with a standard deviation of the pressure greater than 10 bar are, in most cases, caused by air bubbles trapped in one of the piston chambers. To clear trapped air bubbles, it is recommended that you run the MS pump at about 500 $\mu\text{L}/\text{min}$ with methanol or acetonitrile for about 15 minutes. This flow rate is much more effective than purging at higher flow rates with methanol / buffer mixtures.

Strong pump pulsations can also be caused by insufficient flow to the pump. If the above procedure does not reduce the pulsation, replace the solvent reservoir filters and thoroughly purge air from the system.

If it is determined that the cause is not related to trapped air, the cause might be related to one or more damaged piston seals. Replace both piston seals as described in [Chapter 4, "Maintenance."](#)

If the problem still persists after the seals have been changed, the problem is likely due to a contaminated check valve. For low flow ($< 50 \mu\text{L}/\text{min}$) applications, cleaning a check valve is normally not effective. Most likely one or both check valve assemblies needs to be removed and replaced.

Pressure Pulsations with SD between 3 and 10 Bar

Pressure pulsations with a standard deviation of 3 to 10 bar are of various origins. You can continue working, but you should try to find the reason for the pulsation as soon as possible.

Baseline Noise at Low Pressures (under 40 Bar)

HPLC pumps are optimized to deliver a constant flow with a minimum of pressure fluctuations when they are operated in the range of 40 to 400 bar (580 to 5800 psi). The MS pump provides stable flow over the entire pressure range used by HPLC.

Under extreme conditions (flow rates of less than 100 $\mu\text{L}/\text{min}$) and with a low backpressure (less than 40 bar), you might observe a slight instability in the detector baseline at high sensitivity.

This problem can usually be fixed by adding a flow restrictor to your system between the Surveyor MS Pump Plus and autosampler to increase the backpressure to a level greater than 40 bar.

5 Troubleshooting

Insufficient Solvent Flow from the Pump Outlet

Some typical applications susceptible to this problem include the following:

- Flow injection analysis (FIA) using an LC/MS performed at 10 to 50 $\mu\text{L}/\text{min}$ without an HPLC column
- Capillary LC at 3 to 5 $\mu\text{L}/\text{min}$ and 15 to 20 bar

UV detector performance is affected at high sensitivity by any variation of backpressure and flow rate. The Surveyor MS Pump Plus should normally be operated with at least 40 bar of backpressure. If necessary, use a flow restrictor to increase the backpressure of your system.

Insufficient Solvent Flow from the Pump Outlet

If there is insufficient flow when the MS pump is running, the problem is usually a result of one of the following:

- [Pump Cavitation](#)
- [Blocked Proportioning Valve](#)
- [Blocked Check Valve](#)

Pump Cavitation

A restriction in flow from the solvent reservoir to the pump can cause the pump to cavitate, which results in erratic mobile phase delivery. In extreme cases, there might be no flow at all. The reason is usually either a blocked solvent reservoir filter or a restricted solvent inlet line. A possible but much less probable cause could be a restriction inside of the vacuum degasser.

If replacing the solvent reservoir filter does not solve the problem, do the following:

1. Disconnect the solvent reservoir A line from the inlet port of the degasser and verify unrestricted flow of solvent through the tubing.
2. If the solvent flow is normal, reconnect the line to the inlet port of the degasser and disconnect the corresponding degasser outlet line.
3. Check for unrestricted flow of solvent through the degassing chamber. The flow out of the degasser is normally slower than that going to the degasser.
4. Repeat the procedure for solvents B, C, and D.

Suspect a problem with the degasser only if there is virtually no flow coming out of the degasser.

Blocked Proportioning Valve

If flow from the degasser is normal, then the next area to check is the proportioning valve.

❖ To check for blockage in a proportioning valve

1. Set the pump to deliver 100% of solvent A at a flow rate of 1000 $\mu\text{L}/\text{min}$.
2. Measure the actual flow rate.
3. Repeat steps 1 and 2 for each channel.

If one or more of the values are low, clean the LDA by flushing it with the appropriate solvents, as described in “[Cleaning the LDA Components without Disassembling the LDA](#)” on [page 38](#).

Occasionally a valve might freeze closed entirely if it hasn't been used for an extended period of time. To open the valve, insert the valve tool (stainless steel wire) into the valve inlet and push the floating piston away from the valve seat, as described in “[Servicing the Proportioning Valves](#)” on [page 62](#).

Blocked Check Valve

Insufficient flow rate can also be a result of a blockage in one or both check valves. In this case, clean the liquid displacement assembly by flushing it with the appropriate solvents, as described in “[Cleaning the LDA Components without Disassembling the LDA](#)” on [page 38](#).

Irreproducible Retention Times

There are many reasons for retention time problems. Some of the more common causes are:

- An old or damaged column
- Inconsistent stationary phase (Was a new column installed?)
- A validated method lacks robustness
- Changes in mobile phase composition
- Temperature fluctuations
- Pump problems

If you suspect the pump to be responsible, replace the column with a flow restrictor and run a step gradient, as described in “[Testing the Pump Proportioning](#)” on [page 31](#).

Inaccuracy in the steps is most likely caused by contamination of the proportioning valves. Clean the LDA by flushing it with the appropriate solvents, as described in “[Cleaning the LDA Components without Disassembling the LDA](#)” on [page 38](#).

Accessories and Replacement Parts

This chapter contains a list of parts that can be replaced by Thermo Fisher Scientific factory-trained personnel. Accompanying the list of parts are the corresponding part numbers by which they can be ordered from your local Thermo Fisher Scientific service representative.

The replaceable parts are divided into the following sections:

- [Liquid Displacement Assembly \(LDA\) Parts](#)
- [MS Pump Accessory Kit](#)
- [Solvent Interconnect Kit](#)
- [Other Parts and Assemblies](#)

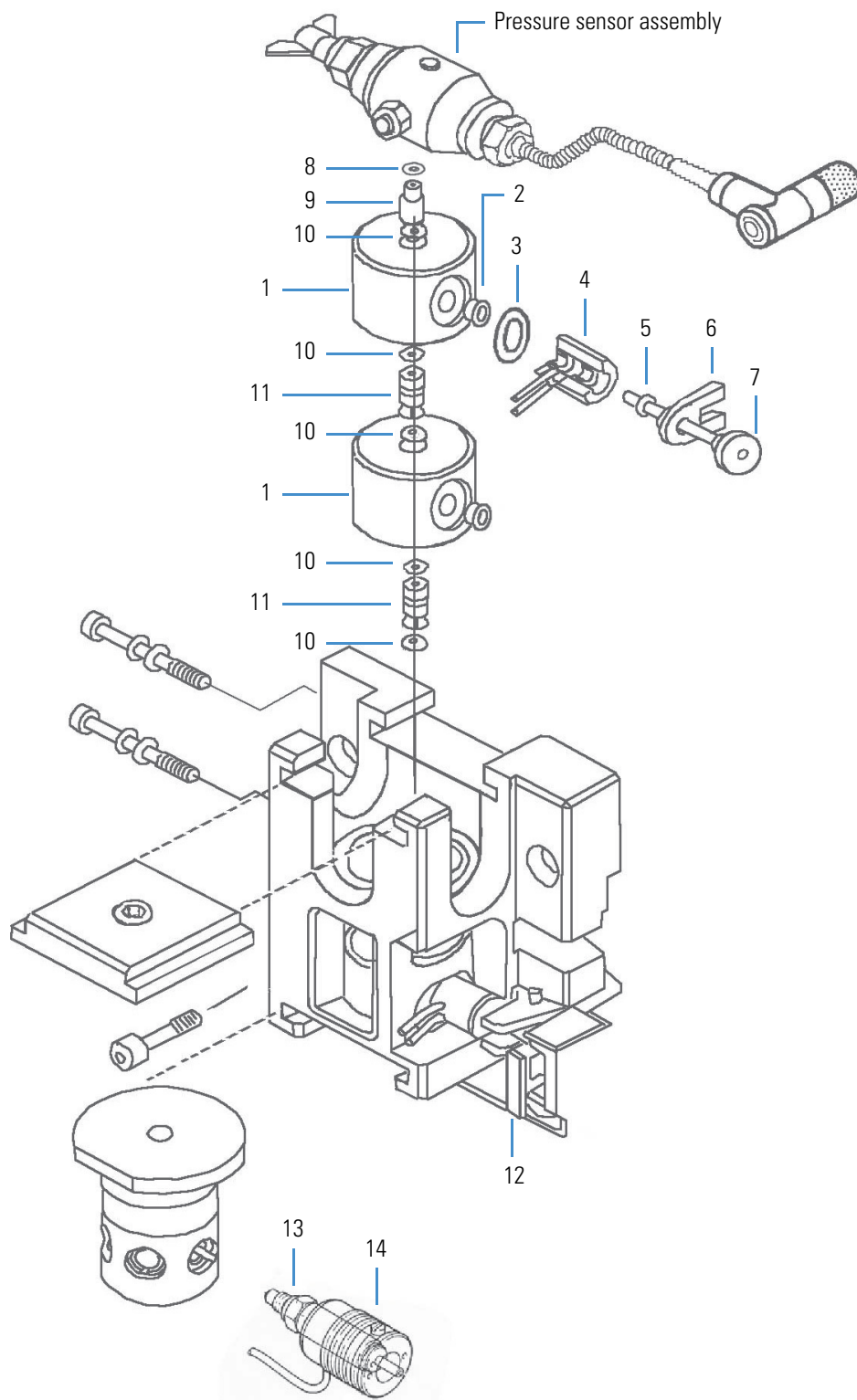
Liquid Displacement Assembly (LDA) Parts

The item numbers in [Table 5](#) refer to [Figure 75](#).

Table 5. LDA Parts

Item No.	Qty	Description	Part Number	Material
1	2	Displacement chamber (piston chamber)	00109-03-00001	Ti
2	2	Primary piston seal	00107-18110	PTFE / GFP / Ti
3	2	Seal Ring, Piston Chamber	60053-01020	PTFE
4	2	Piston guide bushing	00201-11330	SS / Ruby
5	2	Secondary piston seal	00107-18114	PE / SS
6	2	Backing ring	00109-03-00002	SS
7	2	Piston unit	00201-11324	Zirconia
8	1	Outlet cartridge film seal	00107-18112	PEEK
9	1	Outlet cartridge	00109-03-00003	SS
10	5	Film sealing, Face, MS Pump, LDA	00950-01-10013	PEEK
11	2	Check valve cartridge	00110-05110	Ti / Ruby
12	2	Retaining clip	00201-11328	SS
13	4	Proportioning valve	00110--01-00007	Ti / Ruby / Sapphire / Zirconia
14	4	Coil unit for proportioning valve	00201-11344	SS / PEEK
15	4	Sealing ring, proportioning valve	00201-11320	PTFE
Set D	1	Pressure sensor set	00201-11338	PEEK / Ti
	4	Nylon screw, coil	00109-04-00005	Nylon

Figure 75. Liquid displacement assembly



MS Pump Accessory Kit

The following items are contained in the MS pump accessory kit (P/N 72011-62001).

Cable, assembly, RS232, shielded, external PC	72011-63008
Connector, plug, 8 pin, 3.81 mm pitch, Mini-Combicon	00004-02511
Fitting nut / ferrule, Finger-Tight, 1/16 in. OD tube	00101-18088
Fuse, 2 A Slo-Blo (5 × 20 mm) (2 each)	5120-0025
Luer-type adapter	A0796-010
Screwdriver, ball driver, hex, 4 mm	00725-00032
Seal removal tool, P1-4	A3216-010
Seal, H/P, Surveyor MS Pump (2 each)	00107-18110
Syringe, Luer-Lok™, 10 mL	3301-0151
Tool, Prop Valve, MS Pump	60053-01017
Tubing, convoluted, pump waste.	F5034-040
Wrench 1/4 in. and 5/16 in.	5401-0400
Wrench, L-hex, 4 mm	00725-00034

Solvent Interconnect Kit

The following items are included in the LC/MS system solvent interconnect kit.

Bottle cap solvent bottles.	A0343-010
Solvent bottle cap adapter	A3191-010
Solvent filter, 10 micron Teflon.	A4258-010
Tubing, 1/8 in. OD, 240-in length	3219-2004
Fitting, Super Flangeless, 1/8 in. tube, PEEK	00101-18225
Ferrule, Super Flangeless, 1/8 in. tube, PEEK	00101-18017
Label set solvent bottles.	A4158-020
Marker LTR “A”- clip on.	608250001
Marker LTR “B”- clip on.	608250002
Marker LTR “C”- clip on	608250003
Marker LTR “D”- clip on	608250004

Other Parts and Assemblies

The following items include miscellaneous parts and assemblies.

Power cord (European)	6003-0330
Power cord (U.S.A.)	6003-0160
Pulse dampening assembly	00201-11340
RS232 cable	72011-63008
Solvent inlet and outlet line tubing (order by length)	3219-2004
Solvent reservoir filter	A4258-010
Vacuum degas assembly.	00108-00454
Kit, Seal, Surveyor MS Pump	60052-62012
Right door assembly, Surveyor MS Pump Plus	60053-60067
Chassis, Surveyor MS Pump	72011-10001
Front panel, Surveyor MS Pump	72011-10002
PCB assembly, degas monitor and serial	72011-61001
Cable assembly, DI/O to main PCB	72011-63001
Cable assembly, serial to main PCB.	72011-63002
Cable assembly, Degas PCB to main PCB	72011-63003
Cable assembly, DC supply to main PCB	72011-63004
Cable assembly, AC power to Rheos and supply	72011-63005
Cable, motor index	00950-01-00001
Kit, maintenance, Surveyor MS Pump	60052-62010

Tip You can order the individual parts in the No-Ox Tubing Kit. When you order the tubing by the foot, you must cut it to the appropriate lengths. Use no-ox tubing to connect the degassing outlet ports to the proportioning valves.

Tubing, no-ox kit	
(31, 22, 18, and 17 cm lengths of low-pressure tubing and fittings) . . .	00950-01-00021
PEEK fitting.	00101-07-00005
Tefzel ferrule.	0101-08-00010
Kel-F insert.	00101-99-00002
Tubing, no-ox (sold by the foot).	00028-01-00033



Index

Numerics

- 5-connector interconnect cable
 - connections 13
 - figure 12
- 7-connector interconnect cable
 - connections 14
 - figure 14

A

- accessories
 - PEEK fittings 19
 - replacement parts 75
 - stainless steel tubing 5, 19
- accessory kit, contents 10
- adapter cable, MS pump 14
- air bubbles
 - effect on pump pulsation 30, 71
 - effect on retention time reproducibility 67
 - removing 71
 - removing from the piston chambers 37
- autosampler
 - configuring the input signal polarities 23
 - connecting to pump 19

B

- back panel
 - clearance requirements 11
 - connections 11
- backing rings, orientation 56
- backpressure, normal operating 72
- baseline noise, caused by vacuum pump cycling 3
- buffered solutions, pumping 36

C

- cables
 - interconnect 12
 - MS pump adapter cable 14
 - packing location 12
 - power 10
 - RS232 serial communication 2, 12
- capillary LC 72
- captive screws, location 42
- channel, degasser volume 3
- check valves
 - blocked 73
 - cleaning 38, 48
 - contamination 30, 69
 - damaging 47
 - description 3
 - flow direction 52
 - installing 52
 - replacement part 76
 - replacing 47
 - seal rings 48
- checklist 10
- cleaning
 - components of liquid displacement assembly 47
 - pistons 48
- column
 - blockage 66
 - not equilibrated 66–67
- compliance
 - FCC iv
 - regulatory iii
 - WEEE v
- compression lid
 - installing 55
 - removing 43
- compression tool 18
- configuring the signal polarities for the autosampler 23

Index: D

connections
 autosampler 20
 back panel 11
 pulse dampener 20
 solvent lines 15
 system synchronization cable 12
contacting us [xvi](#)
creep caused by salt buildup 36
crossover tube
 function 3
 removing 40

D

DB-9 pin connector 12
degasser
 description 3
 specifications 7
delay volume
 determining 33
 pulse dampener 2
 pump 2–3, 5
 specification 7
detector
 baseline instability 71
 performance 72
documentation
 additional [xiii](#)
 online [xvi](#)
 survey [xvi](#)
dwell volume
 See delay volume

E

electromagnetic compatibility [iii](#)
electronic pulse damping 29

F

FCC compliance [iv](#)
fittings
 connection to autosampler 19
 Super Flangeless 17
flow direction, of check valves 52
flow injection analysis 72
flow rate
 low 71
 minimum and maximum 7
flow restrictor
 used to increase backpressure 71
 using to increase backpressure 31

fuse
 blown 68
 power entry module 68

G

gradient delay volume
 See delay volume

H

holding hooks 57

I

installation
 accessory kit 10
 checklist 10
 serial communication cable 12
 unpacking 10
interconnect cable 12

L

LEDs, front panel 6
Liquid Displacement Assembly (LDA)
 cleaning 47, 73
 components 3, 39
 compression plate 43
 detaching from pump 41
 disassembling 49
 loosening setscrews 43
 schematic diagram 51
low back pressure 71
low flow rates 71

M

maintenance, records 9
manual bleeder valve 27

N

No-Ox Tubing Replacement Kit 79

O

outlet cartridge, installing 54

P

peak area, reproducibility 67
peak shape, effect of plumbing 67
PEEK fittings 10, 19
piston bore 53

- piston chambers
 - installing 54
 - removing 46
- piston guide bushings
 - flushing 36
 - installing 56
 - replacement part 76
- pistons
 - cause of scratches 36
 - cleaning 48
 - movement 58
 - removing 44
 - replacement part 76
 - retaining clips 61
 - salt buildup 36
 - stroke volume 7
- platform, solvent 17
- plunger seals, leaking 69
- power entry module, attaching power cord 21
- power requirements 7
- pressure
 - operating range 7
 - resolution 7
 - settings range 7
 - transducer 3
- pressure readback 30
- pressure sensor assembly
 - installing 55
 - removing 44
 - replacement part 76
- pressure sensor cable
 - connecting 61
 - disconnecting 40
- primary piston seals
 - orientation of spring 50
 - seating 50
- priming the pulse dampener 27
- proportioning accuracy
 - test failure 32
 - testing 31
- proportioning valves
 - blocked 73
 - cleaning 38
 - components 62
 - connecting wires 41
 - contamination 30, 68, 73
 - freezing 63
 - opening 63
 - problems 68
 - replacement part 76

- pulsation test
 - failure 32
 - monitoring backpressure 29
- pulse dampener
 - description 5
 - outlet 27
- pump
 - cable disconnected 68
 - cavitation 72
 - components 2
 - delay volume 3, 5
 - environmental requirements 11
 - leaks 67
 - motor will not run 68
 - powering on 21
 - specifications 7
- pump program, downloading 6
- pump pulsation
 - abnormal 71
 - causes of excessive 30
 - damaged piston seals 71
 - defined 70
 - dirty check valves 71
 - dirty solvent reservoir filters 71
 - effect of trapped air 30
 - monitoring 29
 - normal 70

Q

- quaternary, functional description of pump 2

R

- rear panel connections
 - See* back panel
- records, maintenance 9
- regulatory compliance iii
- retaining clips
 - figure 61
 - installing 61
- retention time, problems 73
- RS232 serial communication cable 12

S

- safety standards iii
- salt buildup
 - cause 36
 - preventing 36
- schematic diagram of the LDA 51
- seal installation tool 50
- seal removal tool 48

Index: T

- seal rings
 - removing 48
 - reusing 49
- seals
 - orientation of primary piston seal 50
 - plunger 69
 - pushing into the piston chamber 50
 - replacement part 76
 - replacing 32, 49
 - ruining 49
 - soaking in methanol 49
- secondary piston seal, replacement part 76
- serial communication
 - configuring the COM port 21
 - connecting the cable 12
- shipping container, contents 10
- signal polarity settings for the autosampler 23
- sinker frit 63
- solvent
 - outlet tubing 19
 - platform 17
- solvent interconnect kit parts 78
- solvent lines
 - connecting 15
 - non-metallic filters 16
 - replacement part 79
- solvent reservoir filter
 - clogged 63
 - replacement part 79
 - replacing 63, 71
- solvents
 - erratic flow 72
 - filtering 15
 - no flow 72
- specifications, pump 7
- status
 - LEDs 2, 6
 - Status page of the Information view 29
- steel compression ring 18
- step-gradient for testing proportioning accuracy 73
- Super Flangeless fittings 17
- survey link xvi

T

- tools
 - 4 mm balldriver wrench 40
 - 4 mm Hex wrench 10, 40
 - Allen wrench 10
 - seal installation 50
 - seal removal tool 48

- troubleshooting
 - high backpressure 66
 - no flow 66
 - no instrument control 67
 - no peaks 66
 - poor peak shape 67
 - retention time reproducibility 67
 - tips 65
 - unstable baseline 66
- Troubleshooting Guide 68

U

- unpacking your pump 10

V

- vacuum degasser
 - description 3
 - specifications 7
- valves
 - check
 - See check valves
 - manual bleeder 27
 - proportioning
 - See proportioning valves

W

- WEEE compliance v

X

- Xcalibur
 - configuring the Surveyor Autosampler 23
 - Direct Control dialog box 25
 - Status page of the Information view 30