



TSQ Series II

TSQ Altis, TSQ Quantis, and TSQ Fortis

Getting Connected Guide

80111-97046 Revision A • June 2018



© 2018 Thermo Fisher Scientific Inc. All rights reserved.

OptaMax NG and TSQ Fortis are trademarks, and Thermo Scientific, TSQ Altis, TSQ Quantis, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

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

The following are registered trademarks in the United States and possibly other countries: Nalgene is a registered trademark of Nalge Nunc International Corporation. Oerlikon Leybold Vacuum is a registered trademark of OC Oerlikon Corporation AG. Rheodyne is a registered trademark of IDEX Health & Science, LLC. SOGEVAC is a registered trademark of Oerlikon Leybold Vacuum. Swagelok is a registered trademark of the Swagelok Company. Tygon is a registered trademark of the division of Saint-Gobain Performance Plastics Corporation.

Chemyx is a trademark of Chemyx Inc. MX Series II is a trademark of IDEX Health & Science, LLC.

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

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: Rev A, June 2018

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 North American and European regulations. Your system meets the applicable requirements in the electromagnetic compatibility (EMC) and product safety standards described in this section.

Unauthorized changes that you make to your system will void regulatory compliance and may defeat the built-in protections for your instrument. Some examples of unauthorized changes include using replacement parts or adding components, options, or peripherals that Thermo Fisher Scientific has not qualified and authorized. Unauthorized changes can also result in bodily injury and/or damage to your system and laboratory.

Ensure continued compliance with regulatory standards:

- Follow all installation instructions provided in the documentation that comes with your system.
- Order replacement parts (as specified in the instrument manual) and additional components, options, and peripherals directly from Thermo Fisher Scientific or an authorized representative.

Low Voltage Directive 2006/95/EC

This device complies with Low Voltage Directive 2006/95/EC and the harmonized safety standard IEC/EN/CSA/UL 61010, 3rd Edition.

EMC Directive 2004/108/EC

This device was tested by TÜV Rheinland of North America and complies with the following EMC standards:

47 CFR 15, Subpart B, Class A: 2012	EN 61000-3-3: 2013	EN 61000-4-8: 2010
CISPR 11: 2009 + A1	EN 61000-4-2: 2009	EN 61000-4-11: 2004
ICES-003 Issue 5: 2012	EN 61000-4-3: 2006 + A1 + A2	BSMI CNS 13438: 2006
EN 55011: 2009 + A1	EN 61000-4-4: 2004 + A1	VCCI V-3: 2015.04
EN 61326-1: 2013	EN 61000-4-5: 2006	
EN 61000-3-2: 2014	EN 61000-4-6: 2013	

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.

WEEE Directive 2012/19/EU



Thermo Fisher Scientific is registered with B2B Compliance ([B2Bcompliance.org.uk](https://www.b2bcompliance.org.uk)) in the UK and with the European Recycling Platform ([ERP-recycling.org](https://www.eur-recycling.org)) in all other countries of the European Union and in Norway.

If this product is located in Europe and you want to participate in the Thermo Fisher Scientific Business-to-Business (B2B) Recycling Program, send an email request to weee.recycle@thermofisher.com with the following information:

- WEEE product class
- Name of the manufacturer or distributor (where you purchased the product)
- Number of product pieces, and the estimated total weight and volume
- Pick-up address and contact person (include contact information)
- Appropriate pick-up time
- Declaration of decontamination, stating that all hazardous fluids or material have been removed from the product

For additional information about the Restriction on Hazardous Substances (RoHS) Directive for the European Union, search for RoHS on the Thermo Fisher Scientific European language websites.

IMPORTANT This recycling program is **not** for biological hazard products or for products that have been medically contaminated. You must treat these types of products as biohazard waste and dispose of them in accordance with your local regulations.

Contents

	Preface	9
	Accessing Documentation	10
	Special Notices, Symbols, and Cautions	10
	Contacting Us	12
Chapter 1	Vacuum System	13
	Connecting the Forepumps to the Mass Spectrometer	14
	Connecting the Forepumps to the Lab Exhaust System	15
Chapter 2	Gas Connections	17
Chapter 3	Syringe Pump and Modular Valve Assembly	19
Chapter 4	Electrical Connections	21
Chapter 5	External Devices	23
	Contact Closure Signal	23
	Communication Connectors	24
Chapter 6	API Source and Waste Container Connections	27
	Installing the API Source	27
	Connecting the Solvent Waste Container	28
Appendix A	Installation Kits and Power Cords	29
	Syringe Pump and Modular Valve Assembly	29
	Single Mechanical Pump Kit	30
	Dual Mechanical Pumps Kit	31
	MS Setup Kit	32
	Power Cords	33
	Index	35

Contents

Preface

The *TSQ Altis, TSQ Quantis, and TSQ Fortis Getting Connected Guide* is intended for the following Thermo Scientific™ triple-stage quadrupole mass spectrometers (MSs):

- TSQ Altis™ (requires two forepumps)
- TSQ Quantis™ (requires one forepump)
- TSQ Fortis™ (requires one forepump)

Use the information in this guide to install the MS system in your laboratory and connect any external liquid chromatography (LC) devices. External devices include those that are controlled from or are independent of Thermo Scientific mass spectrometry applications, such as the Thermo Xcalibur™ data system.

Also, because this guide uses drawings of various connections and component parts to help illustrate procedures, be sure to start from **1**, no matter where it appears.

IMPORTANT Thermo Fisher Scientific field service engineers access some internal components by opening the top cover. Ensure that the instrument location allows for a vertical clearance of at least 92 cm (36 in.) above the instrument. For additional clearance guidelines, refer to the *TSQ Altis, TSQ Quantis, and TSQ Fortis Preinstallation Guide*.

Contents

- [Accessing Documentation](#)
- [Special Notices, Symbols, and Cautions](#)
- [Contacting Us](#)

❖ To suggest changes to the documentation or to the Help

Complete a brief survey about this document by clicking the button below. Thank you in advance for your help.



Accessing Documentation

The TSQ Altis, TSQ Quantis, and TSQ Fortis MSs include complete documentation. For system requirements, refer to the release notes on the software DVD.

❖ To view the product manuals

- (Windows 7) From the Microsoft™ Windows™ taskbar, choose **Start > All Programs > Thermo Instruments > model x.x**, and then open the applicable PDF file.
- (Windows 10) From the Windows taskbar, choose **Start > All Apps > Thermo Instruments > model x.x**, and then open the applicable PDF file.

❖ To view user documentation from the Thermo Fisher Scientific website

1. Go to thermofisher.com.
2. Point to **Services & Support** and click **Manuals** on the left.
3. In the Refine Your Search box, search by the product name.
4. From the results list, click the title to open the document in your web browser, save it, or print it.

To return to the document list, click the browser **Back** button.






Special Notices, Symbols, and Cautions

Make sure you understand the special notices, symbols, and caution labels in this guide. Most of the special notices and cautions appear in boxes; those pertaining to safety also have corresponding symbols. Some symbols are also marked on the instrument itself and can appear in color or in black and white. For complete definitions, see [Table 1](#).




Table 1. Notices, symbols, labels, and their meanings (Sheet 1 of 2)

Notice, symbol, or label	Meaning
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 product.
Note	Highlights information of general interest.
Tip	Highlights helpful information that can make a task easier.

Table 1. Notices, symbols, labels, and their meanings (Sheet 2 of 2)

Notice, symbol, or label	Meaning
	Caution: To prevent possible harm to the user or damage to the product, follow these instructions.
	Chemical hazard: Observe safe laboratory practices and procedures when handling chemicals. Only work with volatile chemicals under a fume or exhaust hood. 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.
	Heavy object: Never lift or move the instrument by yourself; you can suffer personal injury or damage the instrument.
	Risk of eye injury: Eye injury can occur from splattered chemicals, airborne particles, or sharp objects. Wear safety glasses when handling chemicals or servicing the instrument.
	Trip obstacle: Be aware of cords, hoses, or other objects located on the floor.

Contacting Us

Contact	Email	Telephone	QR Code
U.S. Technical Support	us.techsupport.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752	
U.S. Customer Service and Sales	us.customer-support.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752	
Global support	<ul style="list-style-type: none"> ❖ To find global contact information or customize your request <ol style="list-style-type: none"> 1. Go to thermofisher.com. 2. Click Contact Us, select the country, and then select the type of support you need. 3. At the prompt, type the product name. 4. Use the phone number or complete the online form. ❖ To find product support, knowledge bases, and resources <p>Go to thermofisher.com/us/en/home/technical-resources.</p> ❖ To find product information <p>Go to thermofisher.com/us/en/home/brands/thermo-scientific.</p> 		
<p>Note To provide feedback for this document, go to surveymonkey.com/s/PQM6P62 or send an email message to Technical Publications (techpubs-lcms@thermofisher.com).</p>			

Vacuum System

To connect the forepump or forepumps to the MS and the lab exhaust system, follow these procedures.

Contents

- [Connecting the Forepumps to the Mass Spectrometer](#)
- [Connecting the Forepumps to the Lab Exhaust System](#)



CAUTION In addition to reading the forepump's operating and maintenance instructions, do the following:

- Prevent the forepump from overheating by ensuring that there is sufficient air clearance around the pump.
- Maintain the exhaust pressure from atmospheric pressure minus 15 mbar to 1.15 bar absolute (0.15 bar relative).
- Follow the instructions for adding and changing the oil.

1 Vacuum System

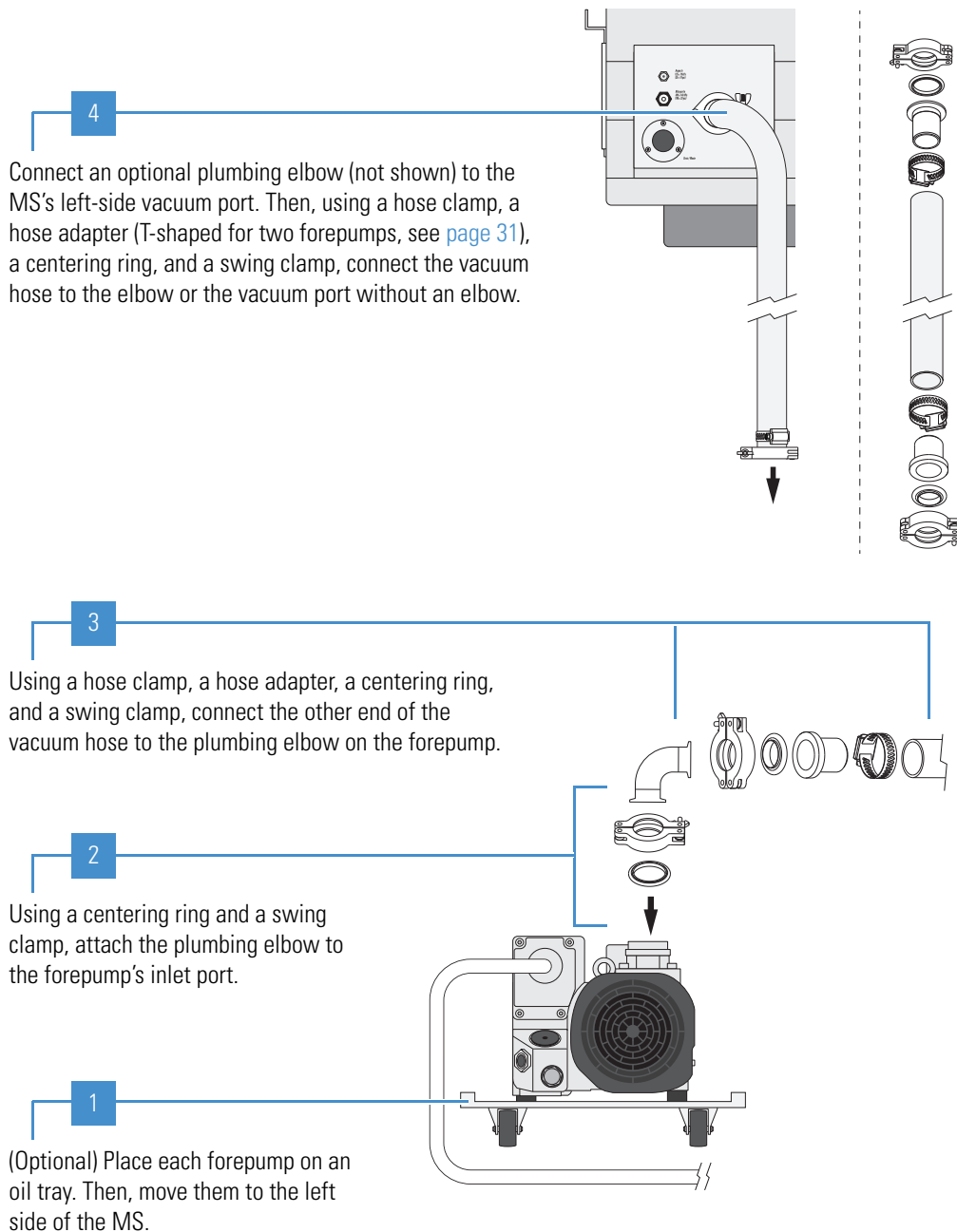
Connecting the Forepumps to the Mass Spectrometer

Connecting the Forepumps to the Mass Spectrometer



IMPORTANT To ensure the best pumping performance, limit the length of the vacuum hose that connects the MS to the forepumps to no more than 2.4 m (8 ft).

❖ To attach the vacuum hose assembly



Connecting the Forepumps to the Lab Exhaust System

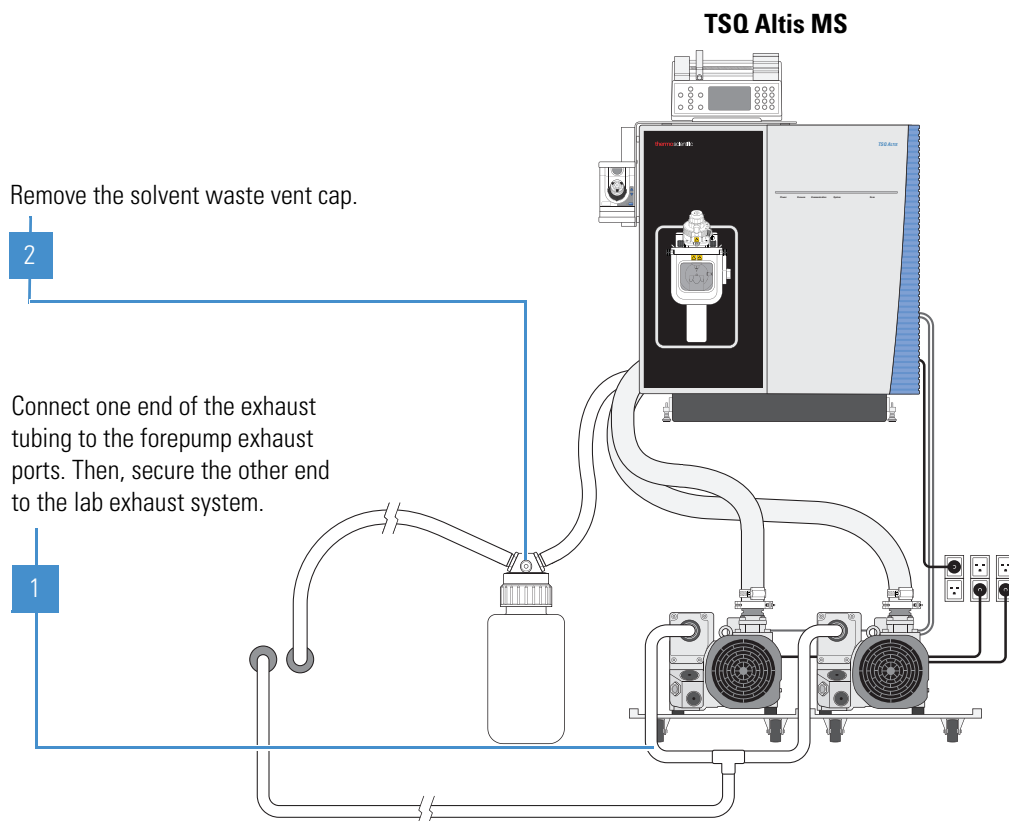
Most atmospheric pressure ionization (API) applications contribute to solvents accumulating in the forepump. Choose an exhaust system that can accommodate the periodic purging of these solvents. The frequency of the purging depends on the throughput of the system—never operate the pump continuously with the gas ballast valve open.



CAUTION

- Run the exhaust hose at floor level for at least 2 m (78.7 in.). This hose acts as a trap for exhaust fumes that would otherwise recondense in the forepump oil.
- The forepump exhaust is a health hazard. Therefore, make sure that the forepump exhaust vents to an appropriate external, active exhaust system.
- Although Thermo Fisher Scientific recommends that you periodically open the gas ballast valve for at least 30 minutes (on the side of the pump) to purge the accumulated solvents, opening the valve might allow a large volume of volatile solvent waste to enter the fume exhaust system.

❖ To connect the forepumps to the exhaust system



1 Vacuum System

Connecting the Forepumps to the Lab Exhaust System

Gas Connections

Before you connect the gas lines to the back of the MS, make sure that the ultra-high-purity (UHP) and high-purity (HP) supply gases are delivered with the necessary pressure and purity as listed in [Table 1](#).

Table 1. Summary of the required gases for the MS

Type	Recommended purity	Required pressure	Typical daily consumption ^a	Function
Argon	UHP (99.995%)—Has less than 1.0 ppm each of water, oxygen, and total hydrocarbons.	135 ±70 kPa (20 ±10 psi)	17 L (0.6 ft ³)	Collision gas
Nitrogen	HP (99%)	690 ±140 kPa (100 ±20 psi)	11 500–26 700 L (406–943 ft ³)	Venting gas; API source auxiliary gas, sheath gas, and sweep gas

^a Approximate value when operating 24 hours and 7 days a week

Note Connections and gas delivery systems might vary. You are responsible for supplying any additional fittings or connections necessary during installation. If your system includes additional devices that require their own gas connections, refer to those instruments' manuals.



CAUTION You must use either copper (provided) or stainless steel for the UHP gas line.



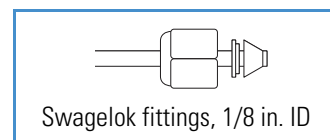
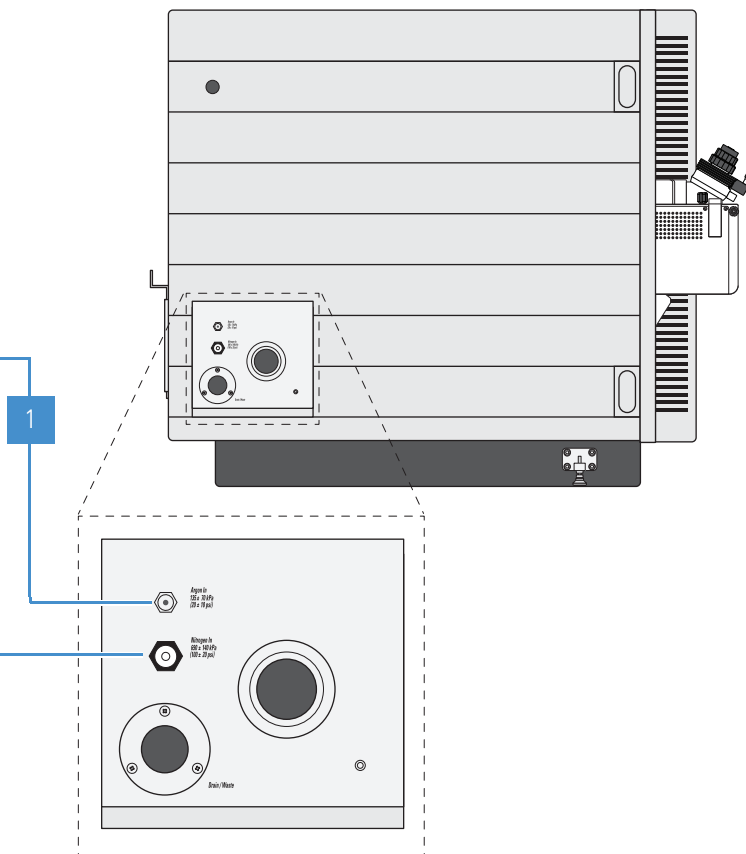
CAUTION After completing these connections, route the tubing so that they are not a trip hazard.

❖ To connect the gas lines to the left side of the mass spectrometer

(UHP argon gas) Cut a length of the copper tubing, attach the 1/8 in. ID Swagelok™-type nut and ferrule set, and then connect it to the **Argon In** inlet until fingertight. On the other end, connect an appropriate fitting that connects to your UHP argon gas supply.

Note Use an open wrench to tighten the fitting on the gas inlet another 1/8- to 1/4-turn.

(HP nitrogen gas) Cut a length of the Teflon™ PFA tubing and push it into the **Nitrogen In** inlet. On the other end, connect an appropriate fitting that connects to your HP nitrogen gas supply.



IMPORTANT

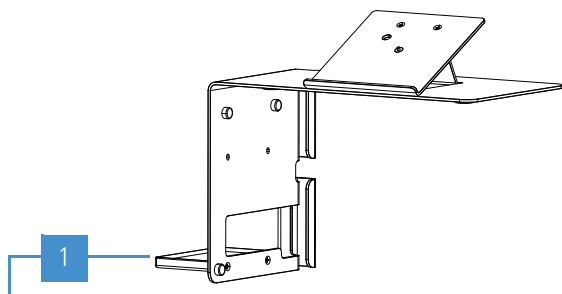
- After you start using the MS, do not shut off the gases. Optimum performance requires a continuous gas flow.
- Do not shut off the UHP nitrogen gas unless you have vented the MS.
- If you intend to use helium for sparging your LC solvents, you must have a second tank and regulator.

Syringe Pump and Modular Valve Assembly

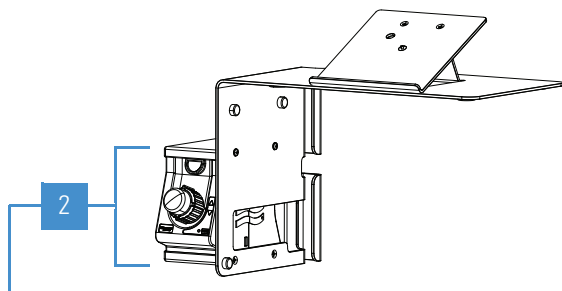
To attach the syringe pump and the modular divert/inject valve, follow this procedure. The only tool required is a #2 Phillips screwdriver.

Note For instructions on how to set up the syringe pump and plumb the modular valve, refer to the *TSQ Altis, TSQ Quantis, and TSQ Fortis Hardware Manual*.

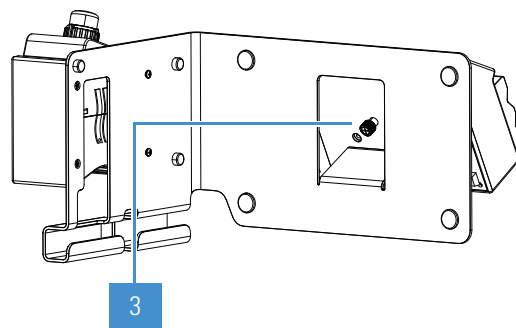
❖ To assemble the syringe pump and modular valve¹



Using the screwdriver and two of the provided screws, attach the bottom shelf to the holder.



Place the valve on the bottom shelf; using the screwdriver and two screws, attach the top shelf.



Center the bottom of the syringe pump on the top angled, support bracket. Then, using your fingers or the screwdriver, tighten the knurled, spring-type screw into the pump.

4

(Not shown) Place the assembly on the top-left surface of the MS and adjust its position as needed.

¹ The holder size might not be as shown.

3 Syringe Pump and Modular Valve Assembly

Electrical Connections

Read the following cautions before you make the electrical connections in the next procedure. If the API source contains the optional APPI lamp, you must provide a USB splitter to accommodate the three USB connections. See also [Chapter 5, “External Devices.”](#)



CAUTION Avoid an electric shock and equipment damage.

- Forepumps—Read the warning on the forepump. When you disconnect the power cord from the forepumps from the electrical outlet do not touch the plug’s pins for at least 5 minutes to allow time for the voltage potential to dissipate.
- Mass spectrometer—Be aware of the following:
 - The AC Output receptacle is not used at this time—do not plug the forepump power cord into this receptacle; see step 5 on [page 10](#).
 - Before you begin the installation, place the Main Power switch in the Off (down) position and the Electronics service switch in the Service Mode (down) position.
- Syringe pump—Make sure that the voltage selector switch is set to the appropriate voltage (110 V or 220 V) for your country or territory.



CAUTION Safety and EMC regulations require the use of Category 5e, shielded Ethernet communication cables, maximum 3 m (10 ft) long.



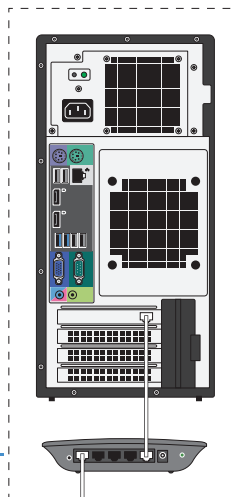
CAUTION After completing these connections, route all cables and power cords so that they are not a trip hazard.

4 Electrical Connections

❖ To connect the MS and data system cables and power cords¹

Connect the **Ethernet cables** from the Ethernet switch to the computer's LC/MS Ethernet network card and the MS.

1

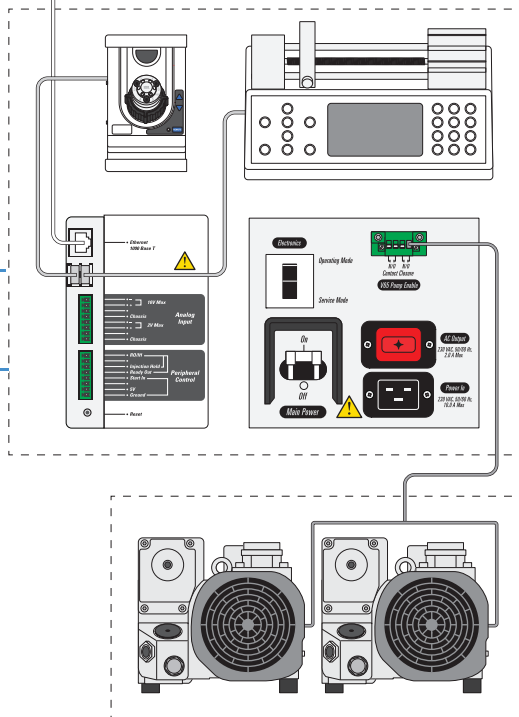


Connect the **USB cables** from the MS to the syringe pump and the modular valve.

2

3

(Not shown) Connect the **contact closure cable** and the **Ready Out cable** (not provided) between the MS and the external device. See Chapter 5, "External Devices."



5

(Not shown) Connect both ends of the **power cords** (for the computer, Ethernet switch, modular valve, syringe pump, MS, and forepumps).

IMPORTANT Plug the power cords for the MS and forepumps into separate, dedicated 230 Vac electrical outlets—that is, circuits A, B, and C for a dual-forepump system.

Note If the modular valve's LED remains off, press its front Remote button to exit local mode.

4

Connect the **relay control cable** between the MS (V65 Pump Enable) and the forepumps.

IMPORTANT Before turning on the MS, make sure to complete all of the system connections described in this guide.

Tip For troubleshooting purposes, you might want to record which devices connect to which ports on the Ethernet switch.

¹ The back of the data system computer might not be as shown.

External Devices

Control of external devices (through a contact closure cable) might or might not be through one of the Thermo Scientific mass spectrometry applications, such as the Xcalibur data system. For instructions on how to set the external device as the start instrument for data acquisition, refer to the *TSQ Altis*, *TSQ Quantis*, and *TSQ Fortis Hardware Manual*.

Contents

- [Contact Closure Signal](#)
- [Communication Connectors](#)

Contact Closure Signal

The MS can start data acquisition upon receiving a contact closure (start) signal from an external LC device, which is typically an autosampler. Thermo Scientific mass spectrometry applications can control external devices (for example, autosamplers, pumps, and detectors) from several manufacturers including Thermo Fisher Scientific, Agilent Technologies, and Waters Corporation.

When an external device is not controlled by a Thermo Scientific mass spectrometry application, such as the Xcalibur data system, you must properly connect it to send its contact closure (start) signal. To assemble the two-wire contact closure cable, use the contact closure mating connector (P/N 00004-21512) supplied in the MS Setup Kit and the contact closure cable provided with the LC device.



CAUTION The external device providing the start signal must have proper earth grounding. Ground loops can cause problems and create a safety hazard. The complementary metal-oxide-semiconductor (CMOS) integrated circuits that are mounted on the internal input/output (I/O) PCB fail if they receive more than 5 Vdc or 5 mA.

IMPORTANT

- Before proceeding, verify that the external device is suitable for use with the MS. Make sure that the signal from the external device meets the requirement stated in the pin 4 (**Start In**) pin-out description (see [Table 1](#)). If it does not meet this requirement, you cannot use it with the MS.
- To help prevent sample waste if the MS becomes unresponsive, connect a cable between the optional LC device and the MS's Ready Out and Injection Hold pins (see [Table 1](#)). Then, configure the Contact Closure settings in the Method Editor. You are responsible for providing this cable.

❖ To connect the contact closure cable

1. Connect the contact closure cable to the **Start In** and Ground pins on the left side of the MS ([Figure 1](#)).
2. Connect the other end of the cable to the external device according to its manual.

Communication Connectors

Use the right-side communication connectors to connect the MS to the Ethernet switch, the syringe pump, the modular valve, and any external device.

[Table 1](#) lists the pin-out descriptions for the connectors shown in [Figure 1](#).

Table 1. Pin-out descriptions for the communication connectors (Sheet 1 of 3)

Pin	Name	Description
–	Reset	Resets the instrument to a power-up state.
<p>Note Use this button only if the instrument does not respond to the control program on the data system computer or if you need to restart the instrument without turning off the electronics service switch.</p>		
Peripheral Control		
1	Ground	Earth ground.
2	5V	Provides a 5 Vdc, 500 mA output (with pin 1).

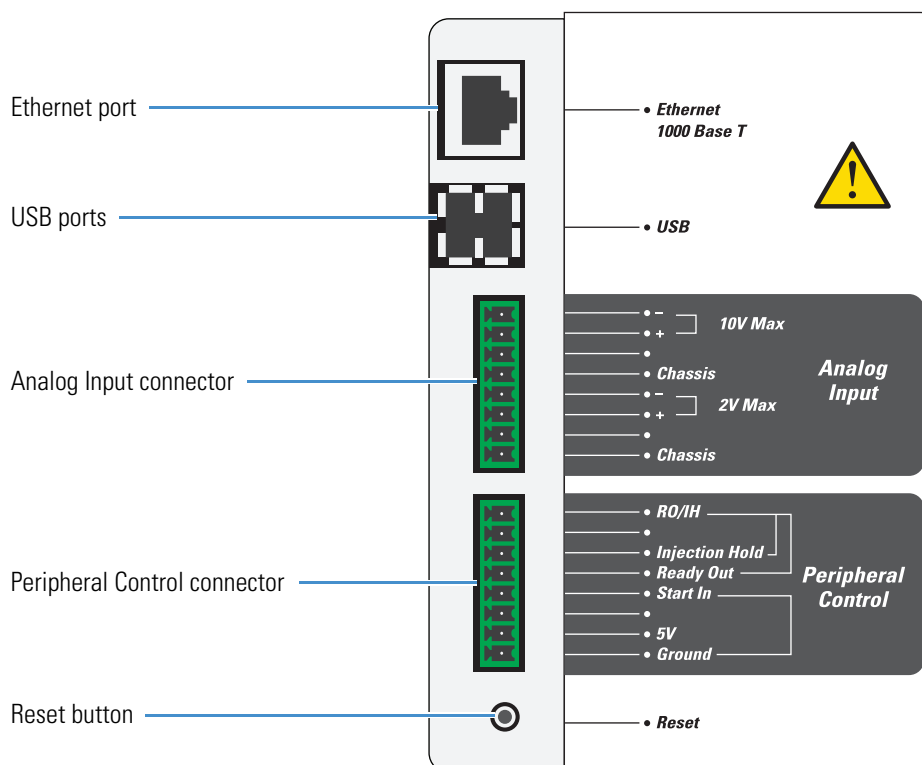
Table 1. Pin-out descriptions for the communication connectors (Sheet 2 of 3)

Pin	Name	Description
4	Start In	<p>Receives the start signal from the contact closure connection of a connected external device.</p> <p>To activate this signal, the external device must pull the signal either low (below 0.75 Vdc) or high (above 2.4 Vdc), depending on the polarity, for at least 100 ms by using a relay, an open-collector driver, or a similar device that connects between pins 4 and 1.</p> <p>Note In the Instrument Configuration window, set the contact closure signal to “High-to-low edge” or “Low-to-high edge,” whichever matches the setting for the connected external device.</p>
5	Ready Out	<p>Provides a relay-driven programmable output signal to the connected external device. The relay opens when a method starts and closes when the method finishes.</p> <p>Output: Maximum 24 Vdc, 3 A</p>
6	Injection Hold	<p>Provides a relay-driven programmable output signal to the connected external device, such as a fraction collector.</p> <p>Output: Maximum 24 Vdc, 3 A</p>
8	RO/IH	Common (return) connection for the Ready Out and Injection Hold pins.
Analog Input		
<p>The two analog channels connect to two separate 12-bit analog-to-digital converters (ADC) for on-demand conversion of the input voltage. The conversion rate depends on the MS rate.</p>		
1	Chassis	Earth ground (for pins 3 and 4).
3, 4	2V Max: + (positive, pin 3) and – (negative, pin 4)	<p>(Channel 2) Provides a connection for an external device, such as an LC instrument.</p> <p>Input: 0–2 Vdc (voltage clamps at 5 Vdc)</p>
5	Chassis	Earth ground (for pins 7 and 8).
7, 8	10V Max: + (positive, pin 7) and – (negative, pin 8)	<p>(Channel 1) Provides a connection for an external device, such as an LC instrument.</p> <p>Input: 0–10 Vdc (voltage clamps at 15 Vdc)</p>

Table 1. Pin-out descriptions for the communication connectors (Sheet 3 of 3)

Pin	Name	Description
Other connectors		
–	USB (2 ports)	Provides a connection for the syringe pump and modular valve.
–	Ethernet 1000 Base T	Provides a connection for the Ethernet switch.

Figure 1. Communication connectors (right side of the MS)



API Source and Waste Container Connections

To install the API source and the solvent waste container to the MS, follow these procedures.

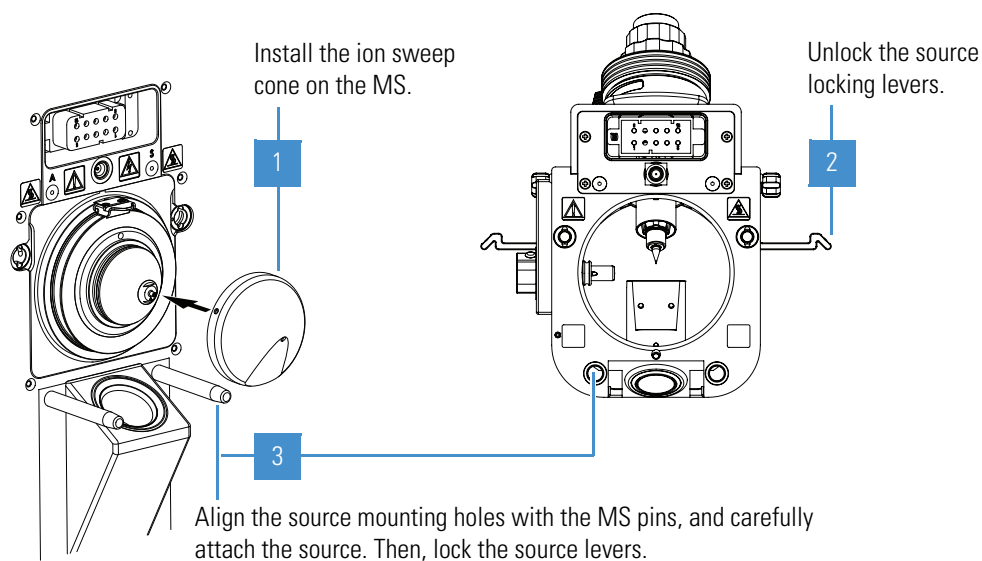
Contents

- Installing the API Source
- Connecting the Solvent Waste Container

Installing the API Source

All the wiring and gas plumbing for the API source are internal. This means you can install or remove the API source or change the ionization mode (H-ESI, APCI, or APPI) by changing the spray insert—all without the use of tools. For information about the spray inserts, optional interface kits, and switching ionization modes, refer to the *OptaMax NG Ion Source User Guide*.

❖ To install the API source onto the MS



Connecting the Solvent Waste Container

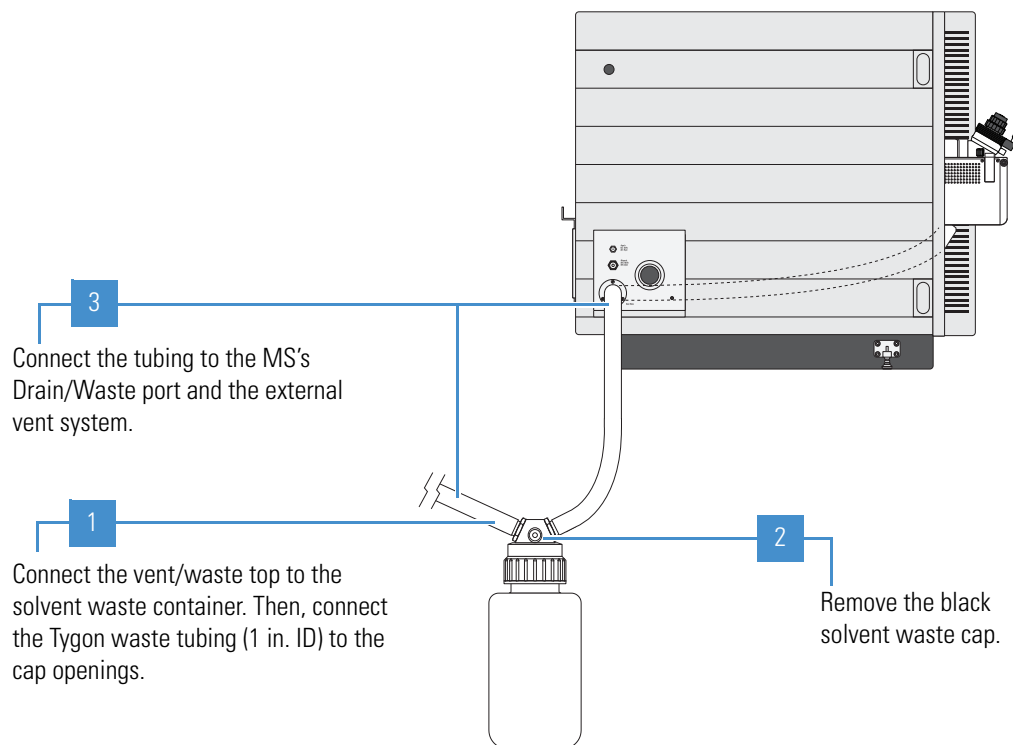
The MS internally routes the solvent waste from the bottom of the API source to the left side drain/waste port. Use the provided Tygon™ tubing to connect the solvent waste container to a fume exhaust system. To prevent solvent waste from backing up into the instrument, make sure that all Tygon tubing is above the level of liquid in the waste container as follows:

- From the MS to the solvent waste container
- From the waste container to the exhaust system

Equip your lab with at least two fume exhaust systems:

- The analyzer optics become contaminated if the drain/waste tubing and the exhaust tubing from the forepump connect to the same fume exhaust system. Route the exhaust tubing from the forepump to a dedicated fume exhaust system.
- Do not vent the Tygon drain tube (or any vent tubing connected to the waste container) to the same fume exhaust system that connects to the forepump. Vent the waste container to a dedicated fume exhaust system. The exhaust system for the API source must accommodate a flow rate of up to 30 L/min (64 ft³/h).

❖ To connect the solvent waste container to the MS



CAUTION After completing these connections, route the tubing so that they are not a trip hazard.

Installation Kits and Power Cords

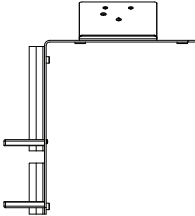
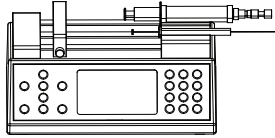
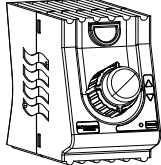
The MS ships with several kits and power cords. This appendix lists the necessary components for the instructions in this guide. For a full list of the kits and their contents, refer to the *TSQ Altis, TSQ Quantis, and TSQ Fortis Hardware Manual*.

Contents

- Syringe Pump and Modular Valve Assembly
- Single Mechanical Pump Kit
- Dual Mechanical Pumps Kit
- MS Setup Kit
- Power Cords

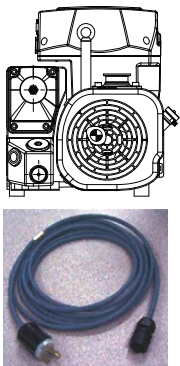

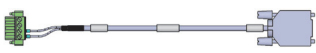
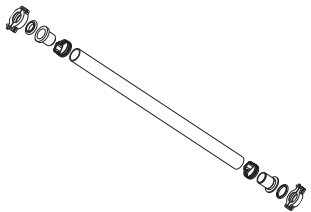
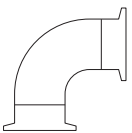

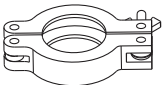

Syringe Pump and Modular Valve Assembly

Table 1. Syringe pump and modular valve assembly

Image	Item	Quantity	Part number
	Holder, L-shaped, syringe pump and modular valve	1	80100-60258
	Chemyx™ Fusion 100T syringe pump with a power cord and USB interface cable	1	00109-99-00045
	Rheodyne™ MX Series II™ modular divert/inject valve (six-port, two-position) with an ac adapter, power cord, USB interface cable, and high-pressure fitting sets	1	00109-99-00046

Single Mechanical Pump Kit

Table 2. Single Mechanical Pump Kit (P/N 80100-62004)

Image	Item	Quantity	Part number
	Forepump, Oerlikon Leybold Vacuum™ SOGEVAC™ SV65BI (rated single-phase 230 Vac, 50/60 Hz)	1	00108-01-00032
	Forepump oil tray, stainless steel, caster wheels	1	00201-99-00549
	Relay control cable, single pump, 2.4 m (8 ft) long (preassembled)	1	80000-63139
	Single pump vacuum hose assembly, KF40 (preassembled; exploded view shown)	1	80000-60229
Elbow connection			
	Elbow, aluminum, NW40, 90 degree	2 ^a	00108-02-00010
	Centering ring with O-ring, nitrile and aluminum, NW40	2 ^a	00108-02-00005
	Swing clamp, aluminum, NW32/40	2 ^a	00108-02-00004
Exhaust waste connection			
	Tubing, Tygon, 3/4 in. (19.1 mm) OD, 0.5 in. (12.7 mm) ID, 3 m (10 ft) long	1	00301-22920

^a Only one piece is required.

Dual Mechanical Pumps Kit

Table 3. Dual Mechanical Pumps Kit (P/N 80100-62013) (Sheet 1 of 2)

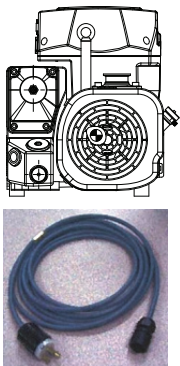

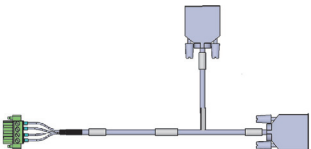
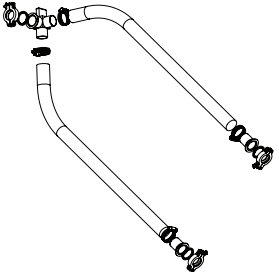
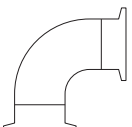

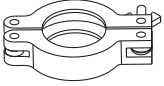


Image	Item	Quantity	Part number
	Forepump, Oerlikon Leybold Vacuum SOGEVAC SV 65 BI	2	00108-01-00032
	Forepump oil tray, stainless steel, caster wheels	2	00201-99-00549
	Dual relay control cable, 2.4 m (8 ft) long (preassembled)	1	80100-63146
	Dual pump vacuum hose assembly (preassembled; exploded view shown)	1	80100-60049
Elbow connection			
	Elbow, aluminum, NW40, 90 degree	3	00108-02-00010
	Centering ring with O-ring, nitrile and aluminum, NW40	3	00108-02-00005
	Swing clamp, aluminum, NW32/40	3	00108-02-00004

Table 3. Dual Mechanical Pumps Kit (P/N 80100-62013) (Sheet 2 of 2)

Image	Item	Quantity	Part number
Exhaust waste connection			
	Fitting Tee, barbed, nylon, for 0.5 in. (12.7 mm) ID tubing	1	00103-01-00012
	Tubing, Tygon, 3/4 in. (19.1 mm) OD, 0.5 in. (12.7 mm) ID, 6 m (20 ft) long	1	00301-22920

MS Setup Kit

Table 4. MS Setup Kit (P/N 80100-62003) (Sheet 1 of 2)




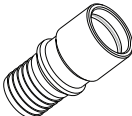

Image	Item	Quantity	Part number
Communications connection			
	Connector plug, MINI-COMBICON™, 8-pin, 26.67 mm (1.05 in.) long, rated 160 V, 8 A (contact closure)	2	00004-21512
–	Connector housing, MINI-COMBICON, 8-position	2	00004-21514
–	Ethernet cables, shielded Category 5e, 2.1 m (7 ft) long	2	00302-99-00036
–	Ethernet power supply (rated 100–240 Vac, 50/60 Hz, 0.6/0.3 A input; 18 W, 12 Vdc, 1.5 A output)	1	00012-01-00039
	Ethernet switch, 5-port Gigabit	1	00825-01-00111
Gas connections			
–	Ferrule, brass, back, 1/4 in. ID	2	00101-04000
–	Ferrule, brass, back, 1/8 in. ID	2	00101-02500
–	Ferrule, brass, front, 1/4 in. ID	2	00101-10000
–	Ferrule, brass, front, 1/8 in. ID	2	00101-08500
–	Swagelok-type nut, brass, 1/4 in. ID	2	00101-12500
–	Swagelok-type nut, brass, 1/8 in. ID	2	00101-15500
–	Tubing, precleaned copper, 1/8 in. OD, 0.030 in. thick, 4.6 m (15 ft) long (for the UHP helium and UHP nitrogen gases)	1 ^a	00301-22701

Table 4. MS Setup Kit (P/N 80100-62003) (Sheet 2 of 2)

Image	Item	Quantity	Part number
–	Tubing, Teflon PFA, 1/4 in. (6.35 mm) OD, 0.062 in. (1.57 mm) thick, 4.6 m (15 ft) long (for the HP nitrogen gas)	1	00101-50100
–	Fittings to connect the copper and PFA tubings to the gas supplies (customer provided)	–	–
Solvent waste connection			
	Container, Nalgene™, 4 L heavy-duty; filling/venting cap	1	80100-20265
	Drain hose adapter with O-ring (not shown)	1	70111-20971
	Tubing, Tygon, 1-3/8 in. OD, 1 in. ID, 3 m (10 ft) long	1	00301-01-00020

^a Your order ships with a second length of copper tubing outside of this kit.

Power Cords

Table 5. Power cord specifications

Device	Cord length	Country or territory ^a	Plug configuration	Plug rating
Mass spectrometer	2.5 m (8 ft)	North America (P/N 96000-98035)	NEMA 6-15	250 Vac, 15 A
		International ^b (P/N 80000-63188)	CEE (3-pole)	250 Vac, 16 A
Forepump	2.5 m (8 ft)	North America (no separate part number)	NEMA 6-15	250 Vac, 15 A
		International (P/N 80000-63186)	CEE (3-pole)	250 Vac, 16 A
Peripheral components	1.8 m (6 ft)	– ^c	–	–

^a Part numbers are provided for Thermo Fisher Scientific field service use only.

^b “International” designation is for countries or territories that do not use the North American plug configuration.

^c Not specified in this guide

A Installation Kits and Power Cords

Power Cords



Index

C

cable connections 22
Communication Connectors 24
communication connectors 26
compliance
 EMC 21
 FCC iii
 regulatory iii
contact closure, start instrument signal 23
contacting us 12
customer-supplied hardware 21, 24

D

directives
 regulatory iii
 WEEE v
documentation, accessing 10

E

electromagnetic compatibility iii
EMC compliance 21
exhaust hose, connecting 15
external devices
 earth ground caution 23
 important notes 24

F

FCC compliance iii
forepump connections
 fume exhaust system cautions 15
 vacuum hose 14
forepumps
 fume exhaust system cautions 28
 gas ballast valve, opening 15
 maintenance instructions 13

G

gases, specifications and consumptions 17

K

kits, installation 29

P

power cords 22, 33
procedures
 API source, installing 27
 contact closure cable, connecting 24
 forepumps, connecting 13
 gases, connecting 18
 modular valve and syringe pump, installing 19
 MS and data system, connecting the cables 22
 solvent waste container, connecting 28

R

regulatory compliance iii
relay switch circuit, external device 25
Reset pin 24

S

safety standards iii
solvent waste system, connecting 28
start signal
 See contact closure, start instrument signal
symbols, meaning 10

T

Thermo Fisher Scientific website, user documents 10
tools 19
troubleshooting, Ethernet switch 22

Index: V

V

valve, LED [22](#)

viewing documents [10](#)

W

WEEE directive [v](#)