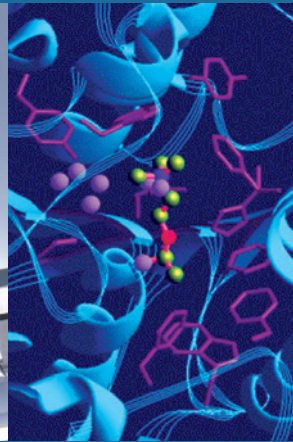
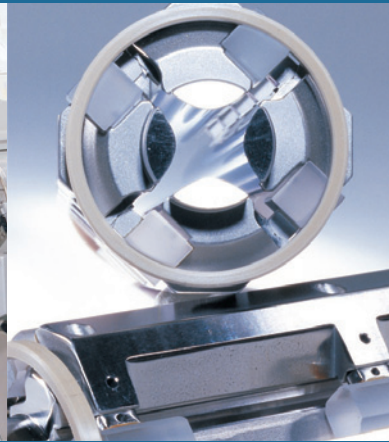


Thermo Fisher Scientific

Orbitrap Velos Pro Hardware Manual

Revision A - 1288290



© 2011 Thermo Fisher Scientific Inc. All rights reserved.

Alconox is a registered trademark of Alconox, Inc. Convectron is a registered trademark of Helix Technology Corporation. Dremel is a registered trademark of the Robert Bosch Tool Corporation. Edwards and the Edwards logo are trademarks of Edwards Limited. HiPace is a trademark of Pfeiffer Vacuum Technology AG. Microsoft and Windows are registered trademarks of Microsoft Corporation. PEEK polymer is a trademark of Victrex plc. PEEKSil is a trademark of SGE International Pty. Ltd. SEQUEST is a registered trademark of University of Washington. Swagelok is a registered trademark of the Crawford Fitting Company. Teflon is a registered trademark of E. I. du Pont de Nemours & Co. Tygon is a registered trademark of Saint-Gobain Performance Plastics Company. Viton is a registered trademark of DuPont Performance Elastomers 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.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

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 released in June 2011.

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

EG-Konformitätserklärung gemäß ISO/IEC 17050-1:2004
EC Declaration of conformity according to ISO/IEC 17050-1:2004

Name des Herstellers: Thermo Fisher Scientific
manufacturers name

Adresse des Herstellers: Hanna-Kunath-Strasse 11
manufacturers address 28199 Bremen
Germany

Der Hersteller erklärt, dass das Produkt

The manufacturer declares that the following product

Name des Produkts: Mass Spectrometer
product name

Modell: ORBITRAP VELOS PRO
model number

Produktoptionen: inkl. / incl. ETD
product options

mit den folgenden EG Richtlinien und harmonisierten Standards übereinstimmt:
is in conformity with the following EC Directives and harmonized standards

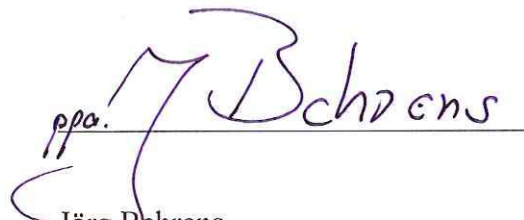
EMV-Richtlinie EMC Directive 2004/108/EG	EN 55011 (11.2007)	EN 61000-4-5 (06.2007)
	EN 61000-3-2 (10.2006)	EN 61000-4-6 (04.2008)
	EN 61000-3-3 (06.2006)	EN 61000-4-11 (02.2005)
	EN 61000-4-2 (12.2001)	EN 61326-1 (10.2006)
	EN 61000-4-3 (12.2006)	+ Corr. (06.2008)
	EN 61000-4-4 (05.2005)	

Niederspannungsrichtlinie EN 61010-1 (08.2002)
Low Voltage Directive
2006/95/EG

Ergänzende Informationen: -
Complementary information

Bremen, Germany, 6. Juni 2011

ThermoFisher
SCIENTIFIC


Jörg Behrens
Technischer Leiter
Director of operations

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.

Velos Pro Mass Spectrometer (April 2011)

EMC Directive 2004/108/EEC

EMC compliance has been evaluated by TUV Rheinland of North America, Inc.

EN 61326-1: 2006	EN 61000-4-3: 2006
EN 55011: 2007, A2: 2007	EN 61000-4-4: 2004
CFR 47, FCC Part 15, Subpart B, Class A: 2009	EN 61000-4-5: 2005
EN 61000-3-2: 2006	EN 61000-4-6: 2007
EN 61000-3-3: 1995, A1: 2001, A2: 2005	EN 61000-4-11: 2004
EN 61000-4-2: 1995, A1: 1999, A2: 2001	

Low Voltage Safety Compliance

This device complies with Low Voltage Directive 2006/95/EEC and harmonized standard EN 61010-1:2001.

Velos Pro/ETD System (April 2011)

EMC Directive 2004/108/EEC

EMC compliance has been evaluated by TUV Rheinland of North America, Inc.

EN 61326-1: 2006	EN 61000-4-3: 2006
EN 55011: 2007, A2: 2007	EN 61000-4-4: 2004
CFR 47, FCC Part 15, Subpart B, Class A: 2009	EN 61000-4-5: 2005
EN 61000-3-2: 2006	EN 61000-4-6: 2007
EN 61000-3-3: 1995, A1: 2001, A2: 2005	EN 61000-4-11: 2004
EN 61000-4-2: 1995, A1: 1999, A2: 2001	

Low Voltage Safety Compliance

This device complies with Low Voltage Directive 2006/95/EEC and harmonized standard EN 61010-1:2001.

FCC Compliance Statement

THIS DEVICE COMPLIES WITH PART 18 OF THE FCC RULES.

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/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on Thermo Fisher Scientific's compliance with these Directives, the recyclers in your country, and information on Thermo Fisher Scientific products which may assist the detection of substances subject to the RoHS Directive are available at www.thermo.com/WEEERoHS.

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 getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliederstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Fisher Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS Anweisung fallen, finden Sie unter www.thermo.com/WEEERoHS.

Conformité DEEE

Ce produit doit être conforme à la directive européenne (2002/96/EC) des Déchets d'Equipements Electriques et Electroniques (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.thermo.com/WEEERoHS.

Read This First

Welcome to the Thermo Scientific Orbitrap Velos Pro system! The Orbitrap Velos Pro is a member of the family of LTQ™ mass spectrometer (MS) hybrid instruments.

All information in this guide concerning the Orbitrap Velos Pro mass spectrometer also applies to the Orbitrap Velos Pro ETD system where the ETD Module is physically coupled to the back of the Orbitrap Velos Pro mass spectrometer.

About This Guide

This *Orbitrap Velos Pro Hardware Manual* contains a description of the modes of operation and principle hardware components of your Orbitrap Velos Pro instrument. In addition, this manual provides step-by-step instructions for cleaning and maintaining your instrument.

Who Uses This Guide

This *Orbitrap Velos Pro Hardware Manual* is intended for all personnel that need a thorough understanding of the instrument (to perform maintenance or troubleshooting, for example). This manual should be kept near the instrument to be available for quick reference.

Scope of This Guide

This manual includes the following chapters:

- **Chapter 1: “Functional Description”** describes the principal components of the Orbitrap Velos Pro mass spectrometer.
- **Chapter 2: “Basic System Operations”** provides procedures for shutting down and starting up the Orbitrap Velos Pro mass spectrometer.
- **Chapter 3: “User Maintenance”** outlines the maintenance procedures that you should perform on a regular basis to maintain optimum instrument performance.
- **Chapter 4: “Replaceable Parts”** lists the replaceable parts for mass spectrometer and data system.

Read This First
About This Guide

- [Appendix A: “Fluoranthene”](#) describes properties of the reagent that is used in the ETD Module portion of the Orbitrap Velos Pro ETD mass spectrometer.

Related Documentation

In addition to this guide, Thermo Fisher Scientific provides the following documents for Orbitrap Velos Pro mass spectrometer and Orbitrap Velos Pro ETD mass spectrometer:

- *LTQ Orbitrap Series Preinstallation Requirements Guide*
- *Orbitrap Velos Pro Getting Started Guide*
- Velos Pro manual set

You can access PDF files of the documents listed above and of this guide from the data system computer. The software also provides Help.

❖ **To view product manuals**

1. From the Microsoft™ Windows™ taskbar, choose **Start > Programs > Thermo Instruments > LTQ > Manuals > *model***.
2. Click the PDF file that you want to view.

Contacting Us

There are several ways to contact Thermo Fisher Scientific.

Assistance

For technical support and ordering information, **visit us on the Web:**

www.thermoscientific.com/ms

Service contact details for customers in Europe are available under:

www.thermoscientific.com/euservicecontact

Customer Information Service

cis.thermo-bremen.com is the Customer Information Service site aimed at providing instant access to

- Latest software updates
- Manuals, application reports, and brochures

Thermo Fisher Scientific recommends that you register with the site as early as possible. To register, visit register.thermo-bremen.com/form/cis and fill in the registration form. Once your registration has been finalized, you will receive confirmation by e-mail.

Changes to the Manual

❖ To suggest changes to this manual

- Please send your comments (in German or English) to:

Editors, Technical Documentation
Thermo Fisher Scientific (Bremen) GmbH
Hanna-Kunath-Str. 11

28199 Bremen

Germany

- Send an e-mail message to the Technical Editor at

documentation.bremen@thermofisher.com

You are encouraged to report errors or omissions in the text or index.
Thank you.

Typographical Conventions

This section describes typographical conventions that have been established for Thermo Fisher Scientific manuals.

Data Input

Throughout this manual, the following conventions indicate data input and output via the computer:

- Messages displayed on the screen are represented by capitalizing the initial letter of each word and by italicizing each word.
- Input that you enter by keyboard is identified by quotation marks: single quotes for single characters, double quotes for strings.
- For brevity, expressions such as “choose **File** > **Directories**” are used rather than “pull down the File menu and choose Directories.”
- Any command enclosed in angle brackets < > represents a single keystroke. For example, “press <F1>” means press the key labeled *F1*.
- Any command that requires pressing two or more keys simultaneously is shown with a plus sign connecting the keys. For example, “press <Shift> + <F1>” means press and hold the <Shift> key and then press the <F1> key.
- Any button that you click on the screen is represented in bold face letters. For example, “click **Close**”.

Topic Headings

The following headings are used to show the organization of topics within a chapter:

Chapter 1 Chapter Name

Second Level Topics

Third Level Topics

Fourth Level Topics

Safety and EMC Information

In accordance with our commitment to customer service and safety, this instrument has satisfied the requirements for the European CE Mark including the Low Voltage Directive.

Designed, manufactured and tested in an ISO9001 registered facility, this instrument has been shipped to you from our manufacturing facility in a safe condition.

This instrument must be used as described in this manual. Any use of this instrument in a manner other than described here may result in instrument damage and/or operator injury.

Notice on Lifting and Handling of Thermo Scientific Instruments

For your safety, and in compliance with international regulations, the physical handling of this Thermo Scientific instrument *requires a team effort* for lifting and/or moving 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: If this instrument is used in a manner not specified by Thermo Fisher Scientific, the protection provided by the instrument could be impaired.

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.

Safety and Special Notices

Make sure you follow the precautionary statements presented in this guide. The safety and other special notices appear different from the main flow of text. Safety and special notices include the following:



Warning Warnings highlight hazards to human beings. Each Warning is accompanied by a Warning symbol. ▲

Caution Cautions highlight information necessary to protect your instrument from damage. ▲

Note Notes highlight information that can affect the quality of your data. In addition, notes often contain information that you might need if you are having trouble. ▲

Identifying Safety Information

This guide contains precautionary statements that can prevent personal injury, instrument damage, and loss of data if properly followed. Warning symbols alert the user to check for hazardous conditions. These appear throughout the manual, where applicable. The most common warning symbols that appear in Thermo Fisher Scientific manuals are shown below.

In addition, every instrument has specific hazards. So, be sure to read and comply with all precautions described in this guide. They will help to ensure the safe and long-term use of your system.



Warning General Hazard. This general symbol indicates that a hazard is present that could result in injuries if it is not avoided. The source of danger is described in the accompanying text. ▲



Warning Electric Shock Hazard. High voltages capable of causing personal injury are used in the instrument. The instrument must be shut down and disconnected from line power before service is performed. Do not operate the instrument with the top cover off. Do not remove protective covers from PCBs. ▲



Warning Burn Hazard. Treat heated zones with respect. Parts of the instrument might be very hot and might cause severe burns if touched. Allow hot components to cool before servicing them. ▲



Warning Corrosive Material. Wear gloves when handling toxic, carcinogenic, mutagenic, or corrosive/irritant chemicals. Use approved containers and procedures for disposal of waste solution. ▲

General Safety Precautions

Observe the following safety precautions when you operate or perform service on your instrument:

- The system should be operated by trained personnel only. Read the manuals before starting the system and make sure that you are familiar to the warnings and safety precautions!
- Accurate results can be obtained only, if the system is in good condition and properly calibrated.

- Service by the customer should be performed by trained qualified personnel only and should be restricted to servicing mechanical parts! Service on electronic parts should be performed by Thermo Fisher Scientific field service engineers only!
- Before plugging in any of the instrument modules or turning on the power, always make sure that the voltage and fuses are set appropriately for your local line voltage.
- Only use fuses of the type and current rating specified. Do not use repaired fuses and do not short-circuit the fuse holder.
- The supplied power cord must be inserted into a power outlet with a protective earth contact (ground). When using an extension cord, make sure that the cord also has an earth contact.
- Do not change the external or internal grounding connections. Tampering with or disconnecting these connections could endanger you and/or damage the system.
- The instrument is properly grounded in accordance with regulations when shipped. You do not need to make any changes to the electrical connections or to the instrument's chassis to ensure safe operation.
- Never run the system without the housing on. Permanent damage can occur. When leaving the system, make sure that all protective covers and doors are properly connected and closed, and that heated areas are separated and marked to protect for unqualified personnel!
- Do not turn the instrument on if you suspect that it has incurred any kind of electrical damage. Instead, disconnect the power cord and contact a Thermo Fisher Scientific field service engineer for a product evaluation. Do not attempt to use the instrument until it has been evaluated. (Electrical damage may have occurred if the system shows visible signs of damage, or has been transported under severe stress.)
- Damage can also result if the instrument is stored for prolonged periods under unfavorable conditions (for example, subjected to heat, water, etc.).
- Always disconnect the power cord before attempting any type of maintenance.
- Capacitors inside the instrument may still be charged even if the instrument is turned off.
- Never try to repair or replace any component of the system that is not described in this manual without the assistance of your Thermo Fisher Scientific field service engineer.

- Do not place any objects upon the instrument—especially not containers with liquids—unless it is requested by the user documentation. Leaking liquids might get into contact with electronic components and cause a short circuit.

Safety Advice for Possible Contamination

Hazardous Material Might Contaminate Certain Parts of Your System During Analysis.

In order to protect our employees, we ask you to adhere to special precautions when returning parts for exchange or repair.

If hazardous materials have contaminated mass spectrometer parts, Thermo Fisher Scientific can only accept these parts for repair if they have been properly decontaminated. Materials that due to their structure and the applied concentration might be toxic or that are reported in publications to be toxic are regarded as hazardous. Materials that will generate synergetic hazardous effects in combination with other present materials are also considered hazardous.

Your signature on the Health and Safety Form confirms that the returned parts have been decontaminated and are free of hazardous materials. Download the form from decon.thermo-bremen.com or order it from the Thermo Fisher Scientific field service engineer.

Parts contaminated by radioisotopes should not be returned to Thermo Fisher Scientific—neither under warranty nor within the exchange part program. If unsure about parts of the system possibly being contaminated by hazardous material, please make sure the Thermo Fisher Scientific field service engineer is informed before the engineer starts working on the system.

Contents

Chapter 1	Functional Description.....	1-1
	General Description	1-2
	Mechanical Characteristics	1-3
	Specifications	1-4
	Control Elements	1-5
	System Status LEDs	1-5
	Control Panels	1-6
	Linear Ion Trap.....	1-12
	Curved Linear Trap	1-13
	Orbitrap Analyzer	1-14
	Extraction of Ion Packets	1-14
	Measuring Principle	1-15
	Ion Detection.....	1-16
	Active Temperature Control	1-17
	HCD Collision Cell.....	1-18
	HCD and ETD	1-18
	ETD System	1-19
	Principle of Operation	1-21
	ETD Module	1-21
	Vacuum System	1-30
	Turbomolecular Pumps.....	1-31
	Forepumps of the Linear Trap	1-33
	Forepump of the ETD Module.....	1-34
	Vacuum System Controls.....	1-35
	Vacuum System Heating during a System Bakeout	1-36
	Gas Supply.....	1-37
	Gas Supply for the Mass Analyzers.....	1-37
	Gas Supply of the Reagent Ion Source	1-40
	Cooling Water Circuit	1-42
	Recirculating Chiller	1-43
	Properties of Cooling Water.....	1-43
	Printed Circuit Boards	1-44
	Linear Ion Trap Electronics.....	1-45
	Electronic Boards on the Right Side of the Instrument	1-46
	Electronic Boards on the Left Side of the Instrument ...	1-59
Chapter 2	Basic System Operations	2-1
	Shutting Down the System in an Emergency	2-2
	Behavior of the System in Case of a Main Failure.....	2-2
	Placing the Instrument in Standby Condition.....	2-4
	Placing the ETD Module in Standby Condition	2-4

	Placing the MS in Standby Condition.....	2-6
	Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely.....	2-7
	Shutting Down the Instrument.....	2-7
	Starting Up the System after a Shutdown.....	2-9
	Starting Up the Instrument.....	2-9
	Setting Up Conditions for Operation.....	2-10
	Starting the ETD Module After a Complete Shutdown	2-11
	Resetting the System	2-12
	Resetting Tune and Calibration Parameters to their Default Values	2-13
	Turning Off the Reagent Ion Source: What to Expect	2-14
Chapter 3	User Maintenance.....	3-1
	General Remarks.....	3-2
	Returning Parts.....	3-3
	Cleaning the Surface of the Instrument.....	3-3
	Maintenance of the Vacuum System	3-4
	Baking Out the System	3-4
	Maintenance of the Forepumps.....	3-5
	Maintenance of the Turbomolecular Pumps	3-11
	Maintenance of the ETD Module.....	3-12
	Handling and Cleaning Reagent Ion Source Parts.....	3-13
	Removing the Access Panels.....	3-18
	Maintenance of the Reagent Ion Source.....	3-20
	Replacing Inlet Valve Components	3-45
	Replacing the Reagent Vials	3-48
	Cleaning the Fan Filters of the ETD Module.....	3-57
	Maintenance of the Cooling Circuit.....	3-58
	Maintenance for the Recirculating Chiller.....	3-58
	Replacing the Water Filter Cartridge.....	3-58
Chapter 4	Replaceable Parts.....	4-1
	Ion Sources	4-2
	Parts for the Basic System.....	4-3
	Parts Lists for the ETD System	4-5
	ETD Reagent Kit.....	4-7
Appendix A	Fluoranthene	A-1
	Glossary	G-1
	Index	I-1

Figures

Orbitrap Velos Pro MS front view	1-2
Schematic of the Orbitrap Velos Pro MS	1-3
Top lid of MS portion opened	1-4
System status LEDs	1-5
Right side of the Orbitrap Velos Pro MS	1-6
Upper control panel	1-7
Power control panel with power control LEDs and switches	1-8
Main power switch	1-9
External connections to the Orbitrap Velos Pro MS	1-10
Layout of the Orbitrap Velos Pro MS, also showing the applied voltages	1-13
Schematic of Orbitrap cell and example of stable ion trajectory	1-14
Principle of electrodynamic squeezing of ions in the Orbitrap analyzer as the field strength is increased	1-15
Approximate shape of ion packets of different m/z after stabilization of voltages	1-16
Schematic of the Orbitrap Velos Pro ETD MS	1-20
Orbitrap Velos Pro ETD MS, rear side	1-22
Rear view of the ETD Module, with major component locations	1-22
ETD Module functional block diagram	1-23
Right side of the Orbitrap Velos Pro ETD MS	1-24
ETD Power Module panel	1-24
Reagent Ion Source dialog box	1-27
Reagent ion source schematics	1-29
Schematic of Orbitrap analyzer vacuum system (CLT compartment and Orbitrap chamber not shown)	1-30
Vacuum components on the left instrument side	1-31
Vacuum components on the right instrument side	1-32
Forepumps cabinet	1-33
Forepump for ETD Module	1-34
Schematic of gas supply for Orbitrap Velos Pro ETD MS	1-37
Proper orientation of the Swagelok-type nut and two-piece ferrule	1-38
Gas regulators	1-39
ETD reagent carrier gas port at the ETD Module	1-40
Schematic of cooling water circuit	1-42
Electronic connections to linear trap	1-45
Electronic boards on the right side of the instrument	1-46
ETD Ion Optic Supply board	1-47
Preamplifier board	1-48
Data Acquisition unit	1-49
Data Acquisition Digital PCI board	1-50
Data Acquisition Analog board	1-51
Instrument Control board	1-52

Power Distribution board	1-54
Power Supply 1 board	1-58
Electronic boards on the left side of the instrument	1-59
Ion Optic Supply board	1-60
Central Electrode Pulser board	1-61
Temperature Controller board	1-62
CLT RF unit (cover removed)	1-64
Central Electrode Power Supply board	1-65
High Voltage Power Supply board (cover removed)	1-67
High Voltage Power Supply board with SPI Bus Termination board	1-68
Main power switch in Off position	2-2
Tune Plus window (Orbitrap Velos Pro ETD), toolbar	2-4
Reagent Ion Source dialog box with Reagent Ion Source On box and Actual condition circled	2-5
Placing the reagent ion source in Standby condition	2-14
Bakeout timer	3-4
Schematic of ETD forepump	3-6
Accessing the ETD Forepump: Removing the panel	3-7
Hooks (left) and top side of detached bottom panel (right)	3-7
Lugs for fixing the bottom panel	3-8
Gas ballast control positions	3-9
Routine maintenance sequence for ETD system	3-12
Rear view of the ETD Module	3-18
Ion source components (left view)	3-21
Tune Plus window (Orbitrap Velos Pro ETD)	3-22
Ion volume tool components	3-23
Guide bar being inserted into guide bar opening	3-23
Guide bar insertion complete	3-24
Rear view of the ETD Module, showing the inlet valve	3-24
Ion volume tool handle in the unlock position	3-25
Ion volume tool guide bar first stop	3-25
Reagent Ion Source dialog box, Open Probe Interlock button	3-26
Instrument Message box: The Ball Valve can now be opened	3-26
Ion volume tool inserted into the inlet valve	3-27
Detail of ion volume tool fully inserted into the inlet valve	3-27
Ion volume tool handle in the locked position	3-28
Ion volume assembly	3-29
Separating ion volume and ion volume holder	3-29
Placing the ion volume on the ion volume tool	3-30
Ion volume tool handle in the unlock position	3-31
Ion volume tool handle in the locked position	3-32
Inlet valve components (ion volume tool not shown)	3-34
Valve shield (1) covering the vacuum manifold probe plate	3-35
Removing the foreline hose from its connection	3-35
Unscrewing the vacuum manifold probe plate	3-36
Removing the vacuum manifold probe plate	3-36
Interior of vacuum manifold	3-37
Removing the ion source assembly from the vacuum manifold	3-38

Ion source assembly 3-39

Ion source assembly exploded view 3-39

Ion source, exploded view 3-41

Ion source lens assembly and ion source 3-43

Filament wire as seen from the bottom of the filament through
the electron lens hole 3-44

Inlet valve components 3-45

Inlet valve seal tool 3-46

Inlet valve seal tool inserted in the inlet valve 3-46

Inlet valve seal on the inlet valve seal tool 3-47

Inlet valve seal disengaged from tool 3-47

Reagent Ion Source dialog box 3-50

ETD Module with back panel removed 3-52

Reagent vials with holders 3-53

ETD Module with vial heater cover removed 3-53

Reagent inlet assembly 3-56

ETD Module, top panel 3-57

Installed water filter 3-59

Removing the filter cartridge 3-60

Filter cartridge with Quick couplers 3-60

ETD Reagent (fluoranthene radical anion) generation from
fluoranthene A-1

Tables

System status LEDs of the Orbitrap Velos Pro MS	1-5
Circuit breakers of the Orbitrap Velos Pro MS	1-7
Typical pressure readings in the ETD Module	1-35
Diagnostic LEDs on the ETD Ion Optic Supply board	1-47
Diagnostic LEDs on the Preamplifier board	1-48
Diagnostic LEDs of the Data Acquisition Digital PCI board	1-50
Diagnostic LEDs of the Data Acquisition Analog board	1-51
Diagnostic LEDs of the Power Supply 2 board	1-52
Diagnostic LEDs of the Instrument Control board	1-53
Software status LEDs of the Instrument Control board	1-53
Status LEDs of the Power Distribution board	1-55
Working modes of the Power Distribution board	1-56
Operating states of the Power Distribution board	1-56
Diagnostic LEDs of the Power Supply 1 board	1-58
Diagnostic LEDs of the Ion Optic Supply board	1-60
Diagnostic LEDs of the Central Electrode Pulser board	1-62
Diagnostic LEDs of the Temperature Controller board	1-63
Diagnostic LEDs of the CLT RF Main board	1-64
Diagnostic LEDs of the Central Electrode Power Supply board	1-66
Diagnostic LEDs of the High Voltage Power Supply board	1-68
User maintenance procedures	3-2
Indications requiring maintenance of the ETD system	3-13

Procedures

To view product manuals.....	iii
To suggest changes to this manual.....	iv
To connect the nitrogen source to the Orbitrap Velos Pro mass spectrometer.....	1-38
To connect the helium source to the Orbitrap Velos Pro mass spectrometer.....	1-38
To place the ETD Module in Standby condition.....	2-4
To place the Orbitrap Velos Pro system in Standby condition.....	2-6
To shut down the instrument completely.....	2-7
To shut down the Orbitrap Velos Pro system.....	2-7
To start up the Orbitrap Velos Pro mass spectrometer.....	2-9
To set up your Orbitrap Velos Pro mass spectrometer for operation.....	2-10
To start up the ETD Module after a complete shutdown.....	2-11
To reset the Orbitrap Velos Pro MS tune and calibration parameters.....	2-13
To perform a system bakeout.....	3-4
To remove the panel.....	3-8
To add oil to the ETD forepump.....	3-8
To purge the rotary-vane pump oil.....	3-9
To change the ETD forepump oil.....	3-10
To clean reagent ion source stainless steel parts.....	3-15
To clean the non-stainless-steel portions of hybrid parts.....	3-17
To remove the ETD main access panel.....	3-18
To remove the ETD side access panel.....	3-19
To clean the ion volume with an inlet valve.....	3-22
To reinsert the ion volume.....	3-29
To clean the ion source lens assembly.....	3-33
To clean the ion source block.....	3-40
To replace the ion source filament.....	3-43
To replace inlet valve components.....	3-45
To place the Orbitrap Velos Pro ETD mass spectrometer in Off Condition and Service mode and to verify that the vials are safe to handle.....	3-49
To install or exchange the reagent vials.....	3-51
To change the reagent ion source flow restrictors.....	3-54
To clean the fan filters of the ETD Module.....	3-57
To replace the water filter cartridge.....	3-59

Chapter 1 Functional Description

This chapter provides an overview of the functional elements of the Orbitrap Velos Pro mass spectrometer. It contains the following topics:

- “General Description” on page 1-2
- “Control Elements” on page 1-5
- “Linear Ion Trap” on page 1-12
- “Curved Linear Trap” on page 1-13
- “Orbitrap Analyzer” on page 1-14
- “ETD System” on page 1-19
- “Vacuum System” on page 1-30
- “Gas Supply” on page 1-37
- “Cooling Water Circuit” on page 1-42
- “Printed Circuit Boards” on page 1-44

General Description

The Orbitrap Velos Pro mass spectrometer is a hybrid mass spectrometer incorporating the Velos Pro™ dual cell linear trap and the Orbitrap™ analyzer. [Figure 1-1](#) shows a front view of the instrument.



Figure 1-1. Orbitrap Velos Pro MS front view

The Orbitrap Velos Pro mass spectrometer consists of four main components (See [Figure 1-2](#) on [page 1-3](#).), which are described in the following topics:

- Dual cell linear ion trap (Thermo Scientific Velos Pro) for sample ionization, selection, fragmentation, and AGC™.
- Intermediate storage device (curved linear trap) that is required for short pulse injection.
- An Orbitrap analyzer for Fourier transformation based analysis.
- Collision cell for performing higher energy CID experiments.

The Orbitrap Velos Pro ETD mass spectrometer has an additional reagent ion source for enabling post-translational modification analyses of peptides by Electron Transfer Dissociation (ETD). See [“ETD System”](#) on [page 1-19](#).

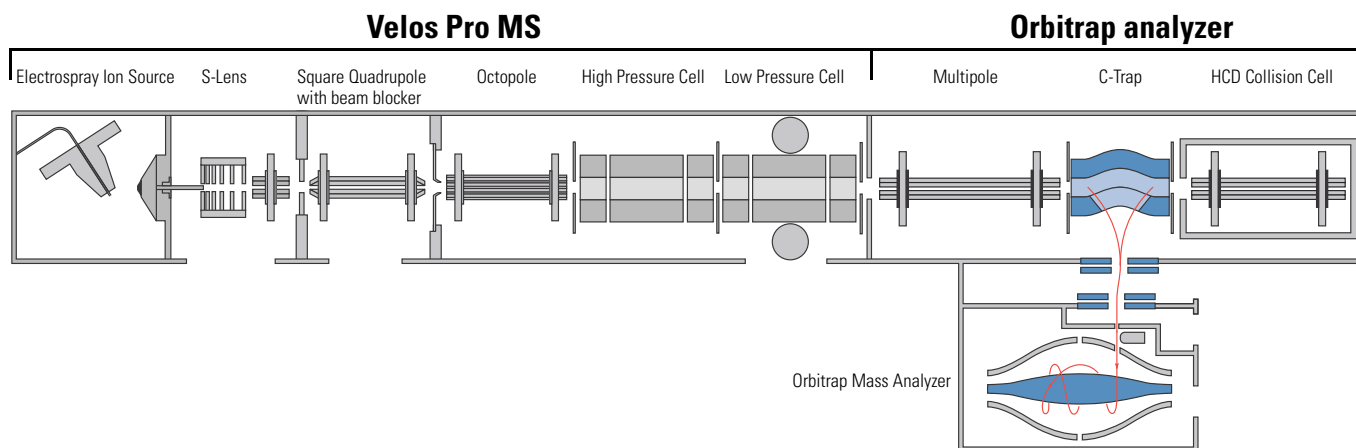


Figure 1-2. Schematic of the Orbitrap Velos Pro MS

Mechanical Characteristics

Wheels at the bottom side of the instrument facilitate positioning the Orbitrap Velos Pro mass spectrometer at the intended place in the laboratory.

The mains inlet as well as a power outlet for peripheral devices are located at the right side of the instrument. The forepumps for the vacuum system of the linear trap and the Orbitrap analyzer are hidden under the linear trap and accessible from the front. The forepump for the ETD Module is accessible after removing the bottom panel of the rear side. The left side panel and the front panel of the MS portion are mounted on hinges and the right side panel is removable. The top lid of the MS portion opens upwards to allow easy access for Thermo Fisher Scientific field service engineers from the top. See [Figure 1-3](#).

In the Orbitrap Velos Pro ETD mass spectrometer, after removing the cables the top lid of the ETD Module is also removable to allow accessing its electronic components.



Figure 1-3. Top lid of MS portion opened

A stand-alone recirculating water chiller is shipped with the instrument. It is connected to the right side of the instrument.

Specifications

The Orbitrap Velos Pro mass spectrometer has the following measuring specifications:

Resolution	60 000 (FWHM) @ m/z 400 with a scan repetition rate of 1 second Minimum resolution 7500, maximum resolution 100 000 @ m/z 400
Cycle Time	1 scan at 60 000 resolution @ m/z 400 per second
Mass Range	m/z 50–2000; m/z 200–4000
Mass Accuracy	<3 ppm RMS for 2 h period with external calibration using defined conditions, <1 ppm RMS with internal calibration
Dynamic Range	>10 000 between mass spectra, >5000 between highest and lowest detectable ion signal in one spectrum
MS/MS	MS/MS and MS^n scan functions

Control Elements

The Orbitrap Velos Pro mass spectrometer is mainly operated from the desktop computer (data system). Some control elements for important system functions are located directly on the instrument. They are described in the following sections.

System Status LEDs

Figure 1-4 shows the system status LEDs at the front of the instrument. Five LEDs indicate the main functions of the system. (See also [Figure 1-5](#) on [page 1-6](#).) The Power LED is directly controlled by the 3 × 230 V input and all other LEDs are controlled by the power distribution board (See [“Power Distribution Board”](#) on [page 1-53](#)). [Table 1-1](#) explains the function of the various LEDs.



Figure 1-4. System status LEDs

The system status LEDs at the front panel of the linear ion trap are described in the *LTQ Series Hardware Manual*.

Table 1-1. System status LEDs of the Orbitrap Velos Pro MS

LED	Status	Information
Power	Green	Main switch on
	Off	Main switch off
Vacuum ^a	Green	Operating vacuum reached
	Yellow	Insufficient vacuum or Vacuum Pumps switch off
Communication	Green	Communication link between instrument and data system established
	Yellow	Communication link starting up or Vacuum Pumps switch off
System ^a	Green	System ready
	Yellow	FT Electronics switch off or Vacuum Pumps switch off
Detect	Blue	Orbitrap analyzer is scanning
	Off	Orbitrap analyzer is not scanning

^aThese LEDs are flashing when a system bakeout is performed. See [“Baking Out the System”](#) on [page 3-4](#).

Control Panels

Figure 1-5 shows the right side of the Orbitrap Velos Pro mass spectrometer. Located here are the control panels, switches, and the ports for the external connections (mains supply, gases, Ethernet communication, and cooling water).

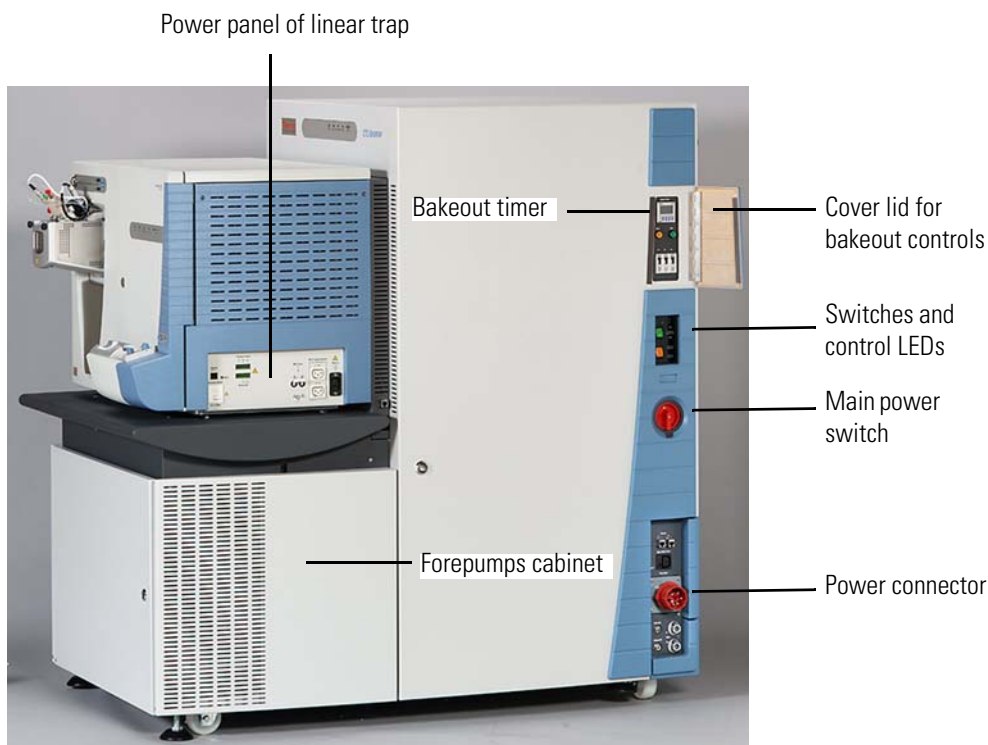


Figure 1-5. Right side of the Orbitrap Velos Pro MS

For more information about the external connections, see “[External Connections](#)” on [page 1-9](#).

Upper Control Panel

The upper instrument control panel comprises the bakeout timer, the bakeout control buttons, and three circuit breakers. To access the upper control panel, swing open the small lid (opens from left to right). See [Figure 1-5](#) and [Figure 1-6](#) on [page 1-7](#).

The timer allows setting the duration for the bakeout of the system. After the duration is set, the bakeout procedure is started by pressing the green button on the right. A running bakeout procedure can be stopped by pressing the orange button on the left side. For instructions about performing a bakeout, see “[Baking Out the System](#)” on [page 3-4](#).

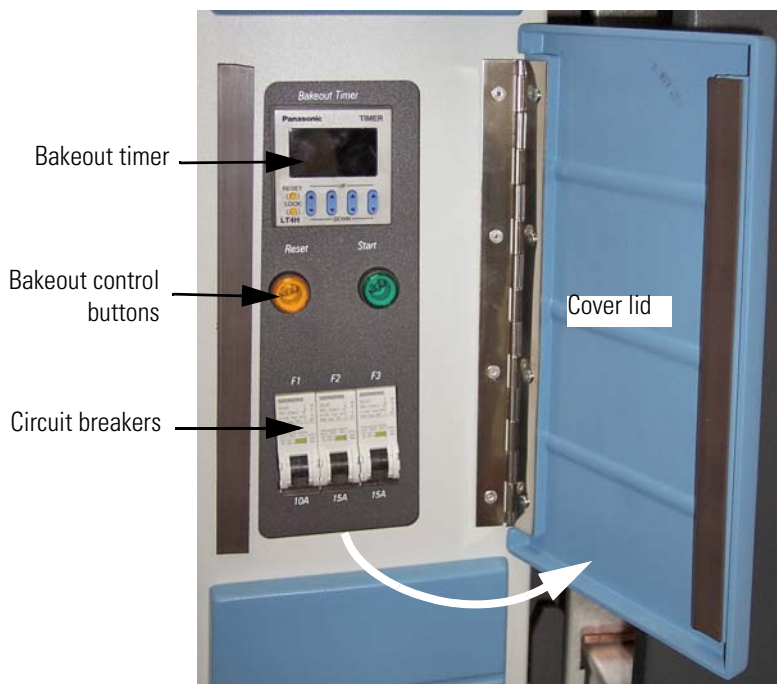


Figure 1-6. Upper control panel

Note The buttons themselves have no indicator function. A running bakeout procedure is indicated by flashing Vacuum and System LEDs at the front side of the instrument. See [Figure 1-4](#) on [page 1-5](#). ▲

Three circuit breakers are located at the bottom of this control panel. [Table 1-2](#) shows the parts of the Orbitrap Velos Pro mass spectrometer that are protected by the respective circuit breaker. The proper function of each circuit breaker is signaled by a dedicated LED in the power control panel (for example, F1 corresponds to L1).

Table 1-2. Circuit breakers of the Orbitrap Velos Pro MS

Circuit breaker	Ampere	LED	Instrument parts
F1	10	L1	Power Distribution
F2	15	L2	Linear ion trap
F3	15	L3	Multiple socket outlets (Peripherals, LC, heater, etc.)

Power Control Panel

In addition to the system status LEDs at the front side (see [Figure 1-4](#) on [page 1-5](#)), the Orbitrap Velos Pro mass spectrometer has three power control LEDs above the Vacuum Pumps switch at the right side. See [Figure 1-7](#). They indicate whether the corresponding circuit breaker is closed and the respective parts of the instrument have power. (See [Table 1-2](#) on [page 1-7](#).)

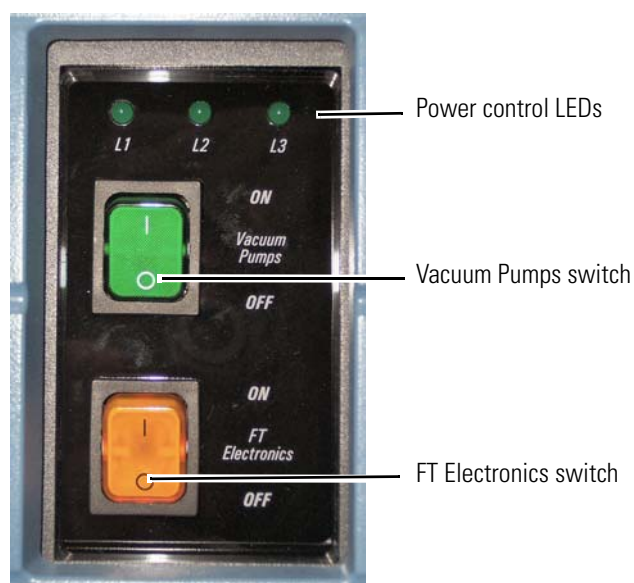


Figure 1-7. Power control panel with power control LEDs and switches

The use of the switches below the power control LEDs changes the working mode of the power distribution. (See [“Working Modes of the Power Distribution”](#) on page 1-48.)

The Vacuum Pumps switch can be set into the positions **ON** or **OFF**. When the switch is in the **OFF** position, everything but the multiple socket outlet is switched off.

The FT Electronics switch can be set into the operating position (**ON**) or into the service position (**OFF**). When the switch is in the service position, all components are switched off with exception of the following:

- Fans
- Heater control
- Power distribution (See [“Power Distribution Board”](#) on page 1-53)
- Pumps (See [“Vacuum System”](#) on page 1-30)
- Temperature controller (See [“Temperature Controller Board”](#) on page 1-62)
- Vacuum control

The linear ion trap also remains on because it has a separate Service switch.

Main Power Switch

The main power switch must be turned 90° clockwise/anti-clockwise to switch on/off the instrument (see [Figure 1-8](#)). Placing the main power switch in the Off position turns off all power to the Orbitrap Velos Pro mass spectrometer, the linear ion trap, and the vacuum pumps. In the Orbitrap Velos Pro ETD mass spectrometer, power to the ETD Module is also turned off.

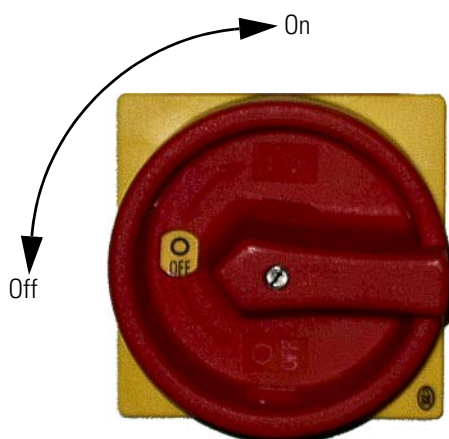


Figure 1-8. Main power switch

Note When the main power switch is in the Off position, you can secure it with a padlock or a cable tie (to prevent unintended re-powering when performing maintenance, for example). ▲

External Connections

[Figure 1-9](#) on [page 1-10](#) shows the lower right side of the instrument with the external connections for mains supply, gases, cooling water, and Ethernet communication.

Located at the top are two ports for Ethernet cables for connecting the Orbitrap Velos Pro mass spectrometer and the linear ion trap via an Ethernet hub with the data system computer.

The power outlet for peripheral devices is located below the Ethernet ports. In the Orbitrap Velos Pro mass spectrometer, the outlet provides the mains supply for the data system. In the Orbitrap Velos Pro ETD mass spectrometer, the outlet provides the mains supply for the ETD Module whereas the data system is connected to a wall outlet.

The power connector for the mains supply is located at the center. The Orbitrap Velos Pro mass spectrometer is designed to operate at a nominal voltage of 230 V AC, 50/60 Hz. Line voltages can vary between a minimum of 207 V AC and a maximum of 253 V AC.

Caution Systems installed in areas with 208 V power experience voltage sags during high use periods that might place the line voltage below the operating parameters discussed in this section. In this case, you must protect your instrument by using a buck/boost transformer to ensure that power is within the specified parameters at all times. ▲

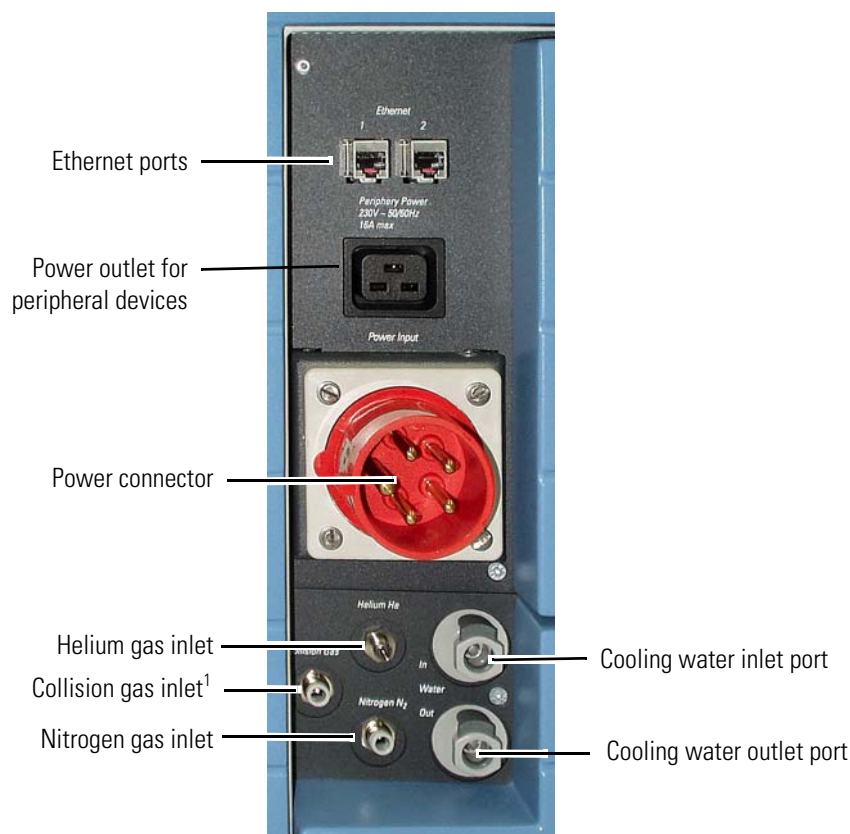


Figure 1-9. External connections to the Orbitrap Velos Pro MS

The cooling water ports are located below the power connector. (See also “Cooling Water Circuit” on page 1-42.)

The port for nitrogen gas allows connecting a Teflon® hose from the gas supply of the laboratory to the instrument. The required gas pressure for nitrogen is 690 ± 140 kPa (6.9 ± 1.4 bar, 100 ± 20 psi). Helium (40 ± 10 psi [275 ± 70 kPa], 99.999% [ultra-high] purity) enters the instrument through a 1/8 inch port. Metal tubing from the helium gas supply must be terminated with 1/8 inch, female, Swagelok-type connectors. See “Gas Supply” on page 1-37 for information about connecting the gas supplies to the instrument.

Caution Do not connect other gases than nitrogen or helium to the Orbitrap Velos Pro mass spectrometer! The maximum pressure for the nitrogen gas inlet is 830 kPa (8.3 bar, 120 psi); the maximum pressure for the helium inlet is 345 kPa (3.45 bar, 50 psi). ▲

¹ The port named Collision Gas is not used in the Orbitrap Velos Pro MS.

In the Orbitrap Velos Pro ETD mass spectrometer, the ETD reagent carrier gas supply of the laboratory is connected via metal tubing to an 1/8 inch inlet port at the rear side of the instrument. Metal tubing from the gas supply must be terminated with 1/8 inch, female, Swagelok-type connectors. The required gas pressure is 690 ± 140 kPa (6.9 ± 1.4 bar, 100 ± 20 psi). See [“Gas Supply of the Reagent Ion Source”](#) on page 1-40.

The exhaust hose from the rotary pumps comes out the back of the instrument, and connects the pumps to the exhaust system in the laboratory.

Linear Ion Trap

The Orbitrap Velos Pro system can utilize a variety of ionization techniques such as ESI, APCI, or APPI. Maintenance of the Ion Max API source, as well as switching between ionization methods, is vent-free. Ions are transferred by octapole and “square” quadrupole lenses into an ion trap that is optimized for axial ion ejection into the curved linear trap. (See [Figure 1-2](#) on [page 1-3](#).)

The linear ion trap is an independent MS detector (Thermo Scientific Velos Pro), which can store, isolate, and fragment ions and then send them either to the Orbitrap analyzer for further analysis or to an SEM detector. The linear ion trap is a unique ion preparation and injection system for Orbitrap MS, because it has greater ion storage capacity than conventional 3D ion trap devices. The linear ion trap is completely described in the *LTQ Series Hardware Manual*.

All the ion handling, selection and excitation capabilities of the ion trap can be used to prepare ions for analysis in the Orbitrap analyzer. These features include storage and ejection of all ions, storage of selected masses or mass ranges, as well as ion isolation. Isolated ions can be excited and then fragmented as necessary for MS/MS and MSⁿ experiments. The patented Automatic Gain Control (AGC) provides extended dynamic range and insures optimized overall performance of the ion trap and Orbitrap MS.

The application of a supplementary RF voltage on the end lenses of the linear trap allows ions of opposite polarity to be trapped in the same space at the same time (charge-sign independent trapping—CSIT). This allows performing ion-ion reactions of previously isolated precursor cations with ETD reagent anions.

The linear ion trap and the transfer chamber are mounted on a table. See [Figure 1-1](#) on [page 1-2](#). The table also serves as a housing for the forepumps. See [Figure 1-25](#) on [page 1-33](#). The Orbitrap Velos Pro mass spectrometer provides power for the linear ion trap. The Orbitrap Velos Pro ETD mass spectrometer also provides the power for the ETD Module.

The linear ion trap is delivered with power connector, gas lines (He, N₂, and collision gas), and vacuum tube lines extending to the ESI source. On the rear side of the Velos Pro ion trap is a flange with an O-ring seal. When the flange is removed, the Orbitrap transfer chamber is mounted to the flange of the linear ion trap. The transfer chamber is held with supports on the table. The components of the ion optics and the Orbitrap analyzer are fixed to the transfer chamber.

Curved Linear Trap

On their way from the linear trap to the Orbitrap analyzer, ions move through the gas-free RF-only octapole into the gas-filled curved linear trap (C-Trap). See [Figure 1-10](#) on [page 1-13](#). Ions entering the C-Trap lose their kinetic energy in collisions with nitrogen bath gas emanating from the HCD cell and get collected near the middle part of the C-Trap. The nitrogen collision gas (bath gas) is used for dissipating the kinetic energy of injected ions and for cooling them down to the axis of the curved linear trap.

Voltages on the end apertures of the curved trap (entrance and exit apertures) are elevated to provide a potential well along its axis. These voltages may be later ramped up to squeeze ions into a shorter thread along this axis. The RF to the C-Trap (“Main RF”) is provided by the CLT RF main board. (See [page 1-63](#).) Entrance and exit DC voltages as well as RF voltages to the octapole are all provided by the ion optic supply board. (See [page 1-60](#).) High voltages to the lenses are provided by the high voltage power supply board. (See [page 1-66](#).)

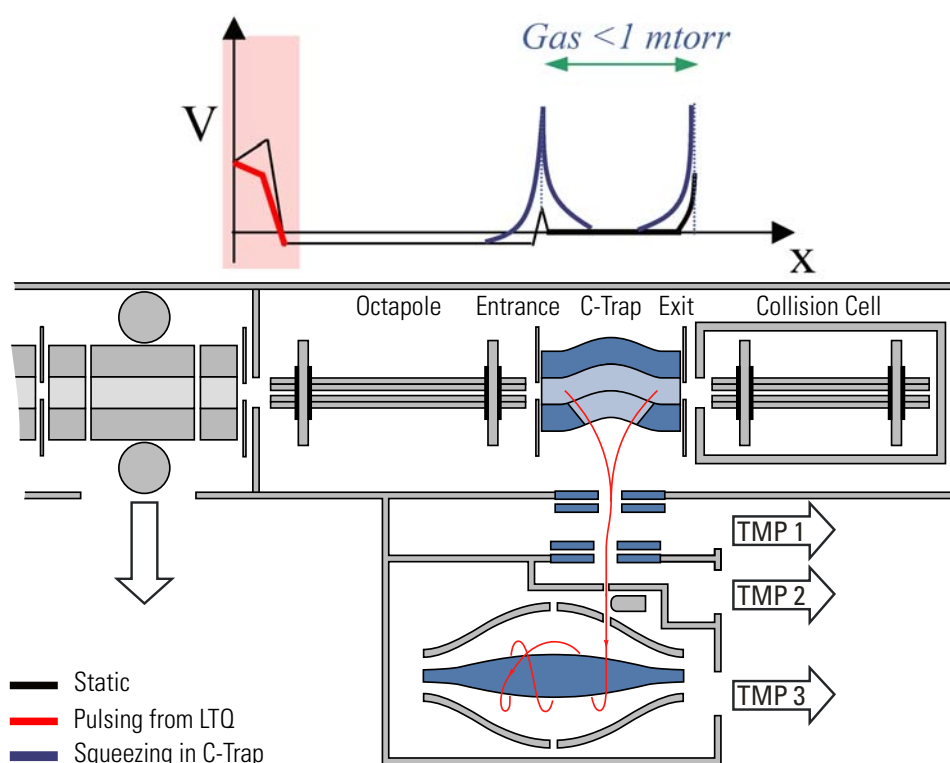


Figure 1-10. Layout of the Orbitrap Velos Pro MS, also showing the applied voltages

Orbitrap Analyzer

The heart of the Orbitrap™ analyzer is an axially-symmetrical mass analyzer. It consists of a spindle-shaped central electrode surrounded by a pair of bell-shaped outer electrodes. See [Figure 1-11](#). The Orbitrap analyzer employs electric fields to capture and confine ions.

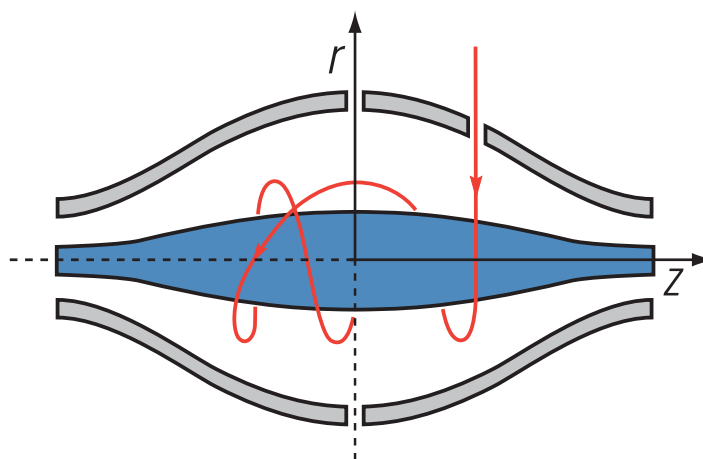


Figure 1-11. Schematic of Orbitrap cell and example of stable ion trajectory

Extraction of Ion Packets

For ion extraction, the RF on the rods of the C-Trap is switched off and extracting voltage pulses are applied to the electrodes, pushing ions orthogonally to the curved axis through a slot in the inner electrode. Because of the initial curvature of the C-Trap and the subsequent lenses, the ion beam converges on the entrance into the Orbitrap analyzer. The lenses that follow the C-Trap (Z-lens) form also differential pumping slots and cause spatial focusing of the ion beam into the entrance of the Orbitrap analyzer. Ions are electrostatically deflected away from the gas jet, thereby eliminating gas carryover into the Orbitrap analyzer.

Owing to the fast pulsing of ions from the C-Trap, ions of each mass-to-charge ratio arrive at the entrance of the Orbitrap analyzer as short packets only a few millimeters long. For each mass-to-charge population, this corresponds to a spread of flight times of only a few hundred nanoseconds for mass-to-charge ratios of a few hundred Daltons/charge. Such durations are considerably shorter than a half-period of axial ion oscillation in the trap. When ions are injected into the Orbitrap analyzer at a position offset from its equator (See [Figure 1-12](#).), these packets start coherent axial oscillations without the need for any additional excitation cycle.

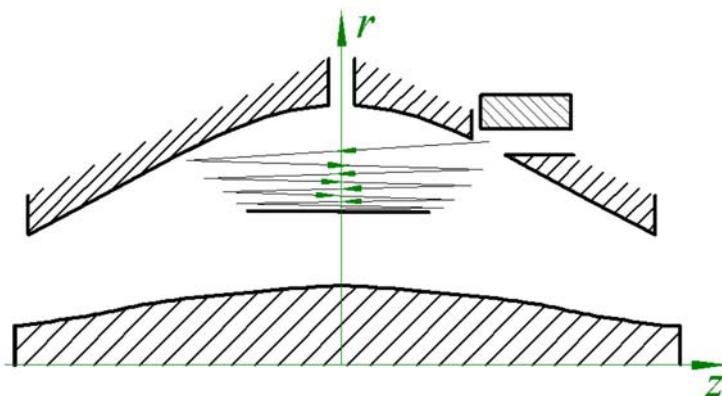


Figure 1-12. Principle of electrodynamic squeezing of ions in the Orbitrap analyzer as the field strength is increased

The evolution of an ion packet during the increase of the electric field is shown schematically on [Figure 1-12](#). When the injected ions approach the opposite electrode for the first time, the increased electric field (owing to the change of the voltage on the central electrode) contracts the radius of the ion cloud by a few percent. The applied voltages are adjusted to prevent collision of the ions with the electrode. A further increase of the field continues to squeeze the trajectory closer to the axis, meanwhile allowing for newly arriving ions (normally, with higher m/z) to enter the C-Trap as well. After ions of all m/z have entered the Orbitrap analyzer and moved far enough from the outer electrodes, the voltage on the central electrode is kept constant and image current detection takes place.

Measuring Principle

In the mass analyzer shown in [Figure 1-11](#) on [page 1-14](#), stable ion trajectories combine rotation around an axial central electrode with harmonic oscillations along it. The frequency ω of these harmonic oscillations along the z -axis depends only on the ions' mass-to-charge ratios m/z and the instrumental constant k :

$$\omega = \sqrt{\frac{z}{m} \times k}$$

Two split halves of the outer electrode of the Orbitrap analyzer detect the image current produced by the oscillating ions. By Fast Fourier Transformation (FFT) of the amplified image current, the instrument obtains the frequencies of these axial oscillations and therefore the mass-to-charge ratios of the ions.

Ion Detection

During ion detection, both the central electrode and deflector are maintained at very stable voltages so that no mass drift can take place. The outer electrode is split in half at $z=0$, allowing the ion image current in the axial direction to be collected. The image current on each of half of the outer electrode is differentially amplified and then undergoes analog-to-digital conversion before processing using the fast Fourier transform algorithm.

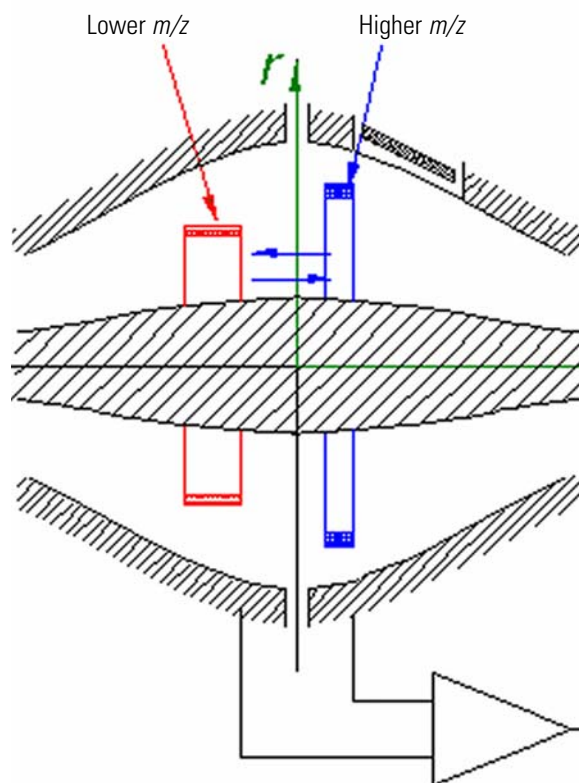


Figure 1-13. Approximate shape of ion packets of different m/z after stabilization of voltages

As mentioned above, stable ion trajectories within the Orbitrap analyzer combine axial oscillations along the z -axis with rotation around the central electrode and vibrations in the radial direction. (See [Figure 1-11](#) on [page 1-14](#).) For any given m/z , only the frequency of axial oscillations is completely independent of initial ion parameters, whereas rotational and radial frequencies exhibit strong dependence on initial radius and energy. Therefore, ions of the same mass-to-charge ratio continue to oscillate along z together, remaining in-phase for many thousands of oscillations.

In contrast to the axial oscillations, the frequencies of radial and rotational motion will vary for ions with slightly different initial parameters. This means that in the radial direction, ions dephase orders of magnitude faster than in the axial direction, and the process occurs in

a period of only 50–100 oscillations. After this, the ion packet of a given m/q assumes the shape of a thin ring, with ions uniformly distributed along its circumference. (See [Figure 1-13](#).) Because of this angular and radial smearing, radial and rotational frequencies cannot appear in the measured spectrum. Meanwhile, axial oscillations will persist, with axial thickness of the ion ring remaining small compared with the axial amplitude. Moving from one half outer electrode to the other, this ring will induce opposite currents on these halves, thus creating a signal to be detected by differential amplification.

Active Temperature Control

Active temperature control is achieved by monitoring temperature directly on the Orbitrap analyzer assembly and compensating any changes in ambient temperature by a thermoelectric cooler (Peltier element) on the outside of the UHV chamber. A dedicated temperature controller board is used for this purpose. See [page 1-62](#).

Peltier Cooling

To allow stable operating conditions in the UHV chamber, it can be cooled or heated (outgassing) by means of a Peltier element located on the outside. A second Peltier element is located on the back of the CE power supply board. See [Figure 1-42](#) on [page 1-59](#).

The Peltier cooling is based on the Peltier Effect, which describes the effect by which the passage of an electric current through a junction of two dissimilar materials (thermoelectric materials) causes temperature differential (cooling effect). The voltage drives the heat to flow from one side of the Peltier element to the other side, resulting in cooling effects on one side and heating effects on the other side.

To remove the heat from the hot side of the Peltier elements, they are connected to the cooling water circuit of the Orbitrap Velos Pro mass spectrometer. See [“Cooling Water Circuit”](#) on [page 1-42](#) for further information.

HCD Collision Cell

The HCD collision cell consists of a straight multipole mounted inside a metal tube, which is connected in direct line-of-sight to the C-Trap. It is supplied with a collision gas to provide increased gas pressure inside the multipole. See “Gas Supply” on [page 1-37](#) for details. The ETD Ion Optic Supply board provides the voltages for the HCD collision cell. (See [page 1-46](#).)

For HCD (Higher Energy Collisional Dissociation), ions are passed through the C-Trap into the HCD collision cell. The offset between the C-Trap and HCD is used to accelerate the parent ions into the gas-filled collision cell. A potential gradient is applied to the collision cell to provide fast extraction of ions, such that it transmits ions at a reliable rate.

The fragment spectra generated in the HCD collision cell and detected in the Orbitrap analyzer show a fragmentation pattern comparable to the pattern of typical triple quadrupole spectra. See the *Orbitrap Velos Pro Getting Started* manual for more information.

HCD and ETD

In the Orbitrap Velos Pro ETD mass spectrometer, ETD reagent anions can efficiently pass through the high pressure region of the HCD cell. This is an important prerequisite to allow for a fast switching (that is, scan to scan) between HCD and ETD fragmentation, thus making comparative measurements possible. When compared with the standard Orbitrap Velos Pro mass spectrometer, HCD performance is not in any way compromised by the addition of the ETD Module.

ETD System

In the Orbitrap Velos Pro ETD mass spectrometer, an ETD Module is physically coupled to the back of the Orbitrap Velos Pro mass spectrometer. A quadrupole mass filter replaces the octapole of the Orbitrap Velos Pro MS. See [Figure 1-2](#) on [page 1-3](#). The linear trap provides the voltages for the quadrupole mass filter. A tube, which contains the transfer multipole, connects the HCD housing to the ETD Module. See [Figure 1-22](#) on [page 1-30](#). The ETD Ion Optic Supply board is mounted on top of the data acquisition unit on the right side of the instrument. See [Figure 1-34](#) on [page 1-47](#).

Protein or peptide analyte ions may also be fragmented in the linear trap by negatively charged reagent ions (fluoranthene radical anions) from the reagent ion source (ETD Module). These negatively charged ions deliver electrons to protein or peptide analyte ions and cause them to fragment at their peptide bonds to produce c and z type fragments (versus the y and b type fragments produced by CID). The resulting analyte fragment ions provide another way of analyzing these molecules as compared to CID and PQD. Electron Transfer Dissociation (ETD) improves the identification of important post-translational modification (PTM) for characterization.

Note Among others, the ETD system is also available as an upgrade on new and existing Velos Pro and Orbitrap Velos Pro systems. ▲

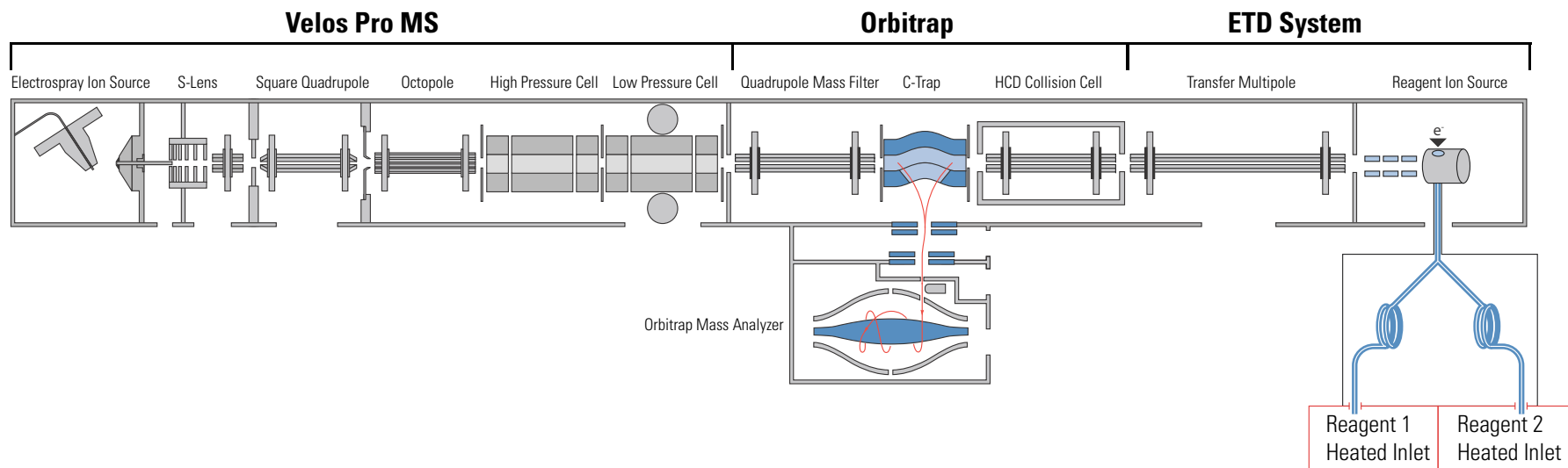


Figure 1-14. Schematic of the Orbitrap Velos Pro ETD MS

Principle of Operation

During a typical ETD MS/MS scan, analyte cations are injected into the linear trap for subsequent precursor cation isolation. Then, ETD reagent anions are generated in the CI ion source and are transferred into the linear trap via RF-only ion guides (transfer multipoles), the gas-filled HCD collision cell, and the C-Trap. (See [Figure 1-14](#) on [page 1-20](#).)

The reagent ions pass a quadrupole mass filter between C-Trap and linear trap. This ion guide works as a low pass mass filter to remove the adduct of the fluoranthene radical and molecular nitrogen at m/z 216. This adduct favors proton transfer reactions instead of electron transfer.

The application of a supplementary RF voltage on the end lenses of the linear trap allows ions of opposite polarity to be trapped in the same space at the same time (charge-sign independent trapping—CSIT).

During ion-ion reactions in the linear trap, electrons are transferred from the reagent anions to the precursor cations. The resulting product ions are mass-to-charge (m/z) analyzed in either the linear trap (if speed and sensitivity are important) or the Orbitrap analyzer (if mass resolution and accuracy are important).

ETD Module

[Figure 1-15](#) on [page 1-22](#) shows the rear side of the ETD Module. It consists of the reagent ion source, ETD Module electronics, ETD Module power supply, ETD Module forepump, and the hardware that connects the ETD Module to the mass detectors.

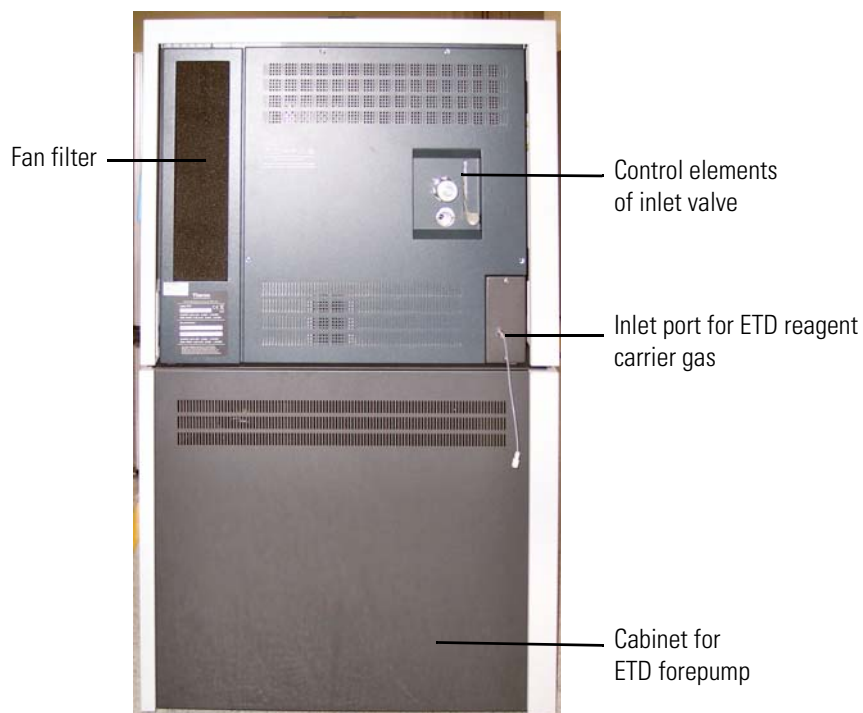
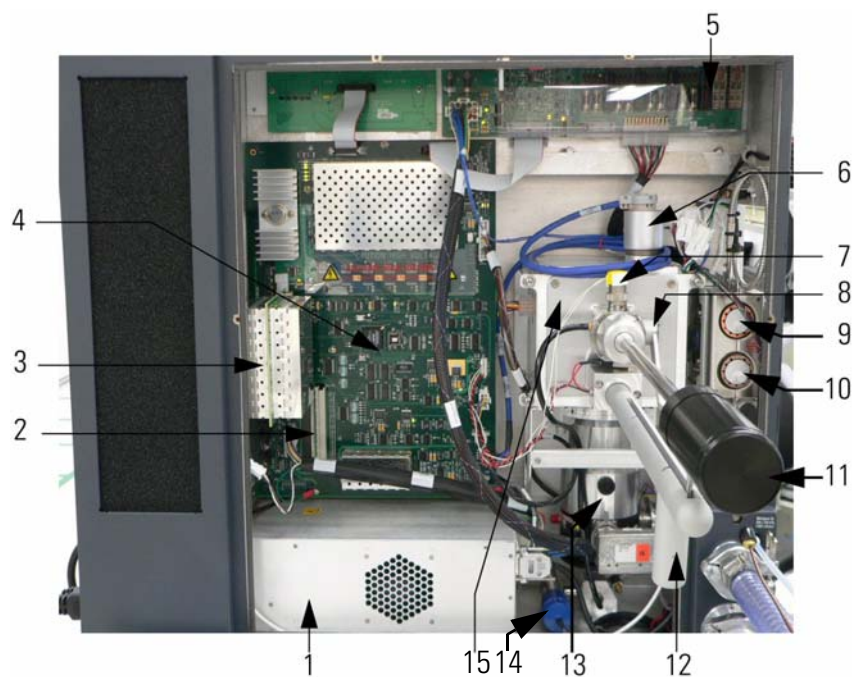


Figure 1-15. Orbitrap Velos Pro ETD MS, rear side



Labeled components: 1=power module, 2=connector to Interface Board (Interface Board is behind the ETD Control PCB, item #4), 3=DC HV Supply PCB, 4=ETD Control PCB, 5=Heater Control PCB, 6=ion gauge, 7=inlet valve solenoid, 8=inlet valve lever in closed (down) position, 9=reagent heater 1, 10=reagent heater 2, 11=ion volume tool handle, 12=guide bar, 13=TMP, 14=Convectron™ gauge, 15=vacuum manifold (contains ion source and ion volume)

Figure 1-16. Rear view of the ETD Module, with major component locations

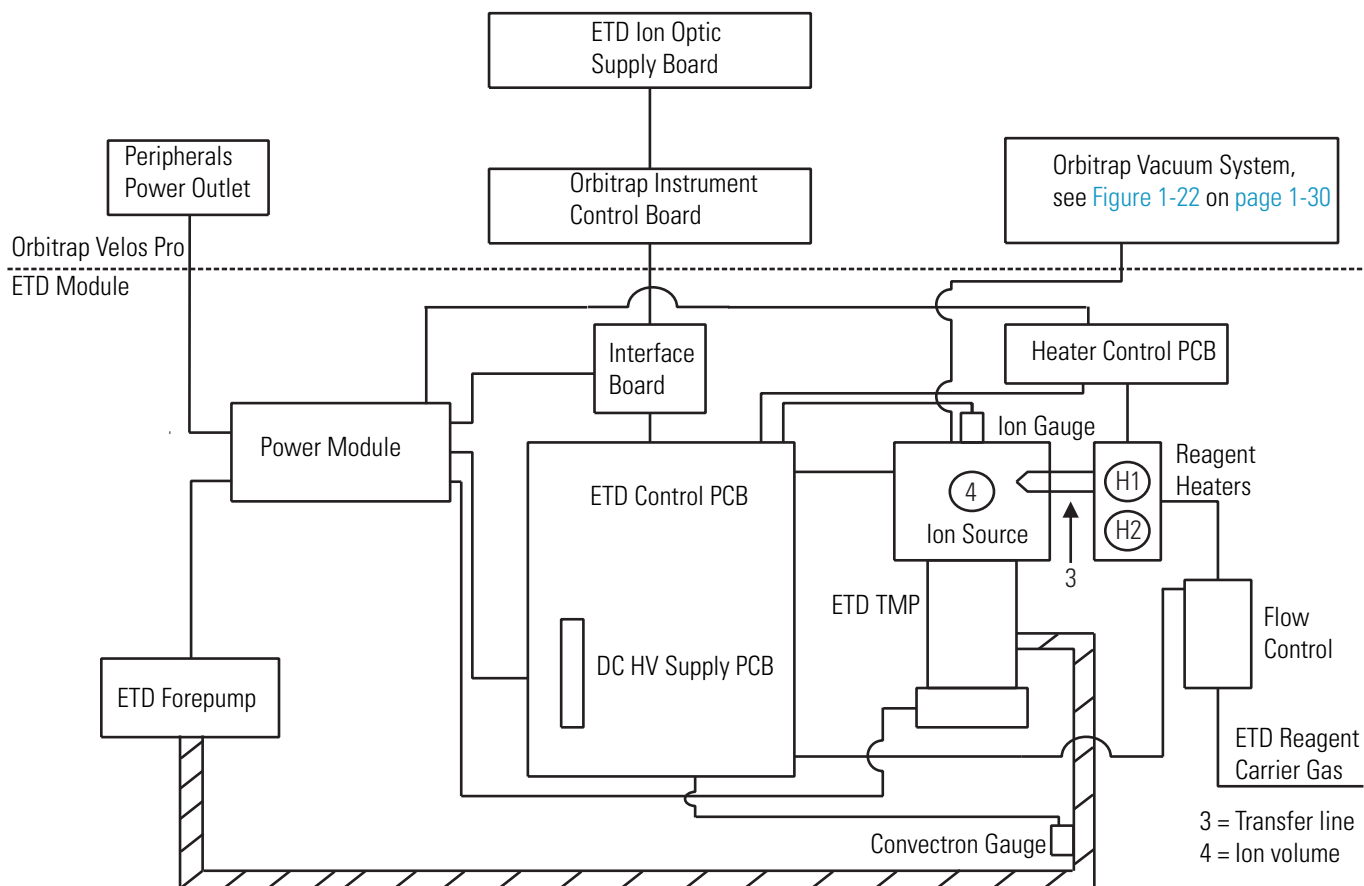


Figure 1-17. ETD Module functional block diagram

The following sections describe the major ETD Module components that are shown in [Figure 1-16](#) on [page 1-22](#) and [Figure 1-17](#).

ETD Power Module

The ETD power module (item #1 in [Figure 1-16](#)) receives 220 V, 10 A, from the peripherals power outlet. See [Figure 1-9](#) on [page 1-10](#). It distributes appropriate voltages and currents to the ETD components. It also contains DC power supplies.

ETD Module Power Panel

The external receptacles and switches for the power module are located on the ETD power module panel at the right side of the ETD Module. See [Figure 1-18](#).



Figure 1-18. Right side of the Orbitrap Velos Pro ETD MS

Figure 1-19 shows a close up picture of the ETD Power Module panel. Power In is connected to the peripherals power outlet of the MS portion. See Figure 1-9 on page 1-10. Forepump is a receptacle to power the ETD forepump (220 V AC, 5 A).

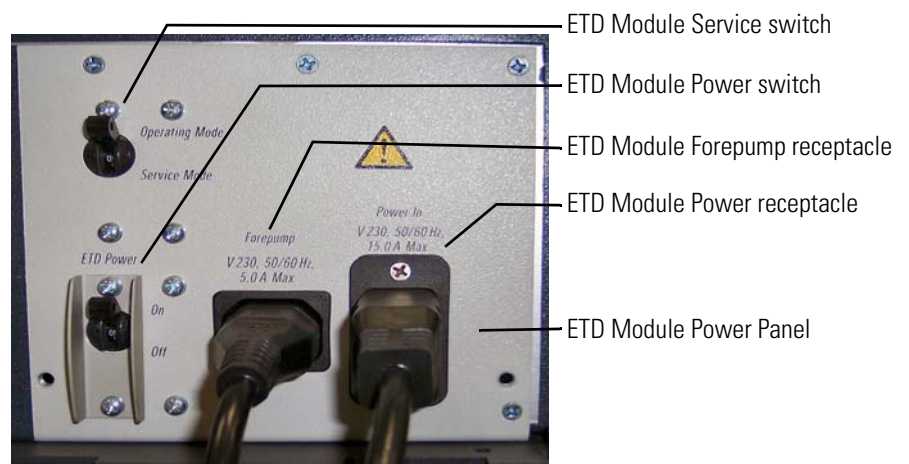


Figure 1-19. ETD Power Module panel

The ETD power module panel contains the main breaker and the service switch for the ETD Module. During normal operation, the ETD Power switch is left On and the service switch is left in the Operating Mode position. As a safety feature, both components of the Orbitrap Velos Pro ETD system (the mass spectrometer and the ETD Module) are shut down with one set of switches, the mass spectrometer switches. When you perform maintenance on components inside the ETD Module as described in “[Maintenance of the ETD Module](#)” on [page 3-12](#), you set the mass spectrometer’s service switch to the Service position. The service switch turns On or Off power to all ETD Module components except turbomolecular pump and forepump.

ETD Module Interface Board

The ETD Module Interface board (item #2 in [Figure 1-16](#) on [page 1-22](#)) provides an electronic interface between the ETD Module and the MS. This board also allows the power to both the MS and the ETD Module to be controlled by the MS power control panel switches:

- The MS Main Power switch controls the power supply to all components in both the MS and the ETD Module.
- The MS FT Electronics switch controls the power supply to all mass spectrometer and ETD Module components except the pumps that are connected to the MS and the ETD Module.

Note The ability to control the power to both components of the Orbitrap Velos Pro ETD mass spectrometer at one point (the power control panel switches of the MS) is a safety feature. ▲

ETD Control PCB

The ETD Control PCB (item #4 in [Figure 1-16](#)) controls most of the ETD Module functions. The ETD Control PCB consists of circuits that control:

- ETD Module operating logic
- Ion source (filament, ion source heater, lenses)
- Heater temperature and readback logic (for reagent heaters, transfer line heater, and the restrictor oven heater)
- Reagent gas flow
- Oven cooling gas control
- Ion gauge
- Convectron™ gauge

The DC HV Supply PCB (item #3 in [Figure 1-16](#) on [page 1-22](#)) is plugged in to the ETD Control PCB.

ETD Heater Control PCB

The ETD Module Heater Control PCB (item #6 in [Figure 1-16](#)) contains the power source and temperature sensing circuitry for the four heaters in the reagent ion source. The heaters are H1 and H2 (the two reagent heaters, [Figure 1-17](#) on [page 1-23](#), and items #9 and #10 in [Figure 1-16](#)), the transfer line heater (#3 in [Figure 1-16](#)), and the restrictor oven heater (not shown in [Figure 1-16](#)). The Heater Control PCB reports temperature information to the heater temperature and readback logic on the ETD Control PCB. The heater temperature and readback logic controls how the Heater Control PCB applies power to the ETD Module heaters.

Reagent Carrier Gas Flow Control for ETD

The ETD Module contains a digital flow control for the chemical ionization (CI) gas/reagent carrier gas provided by the ETD Control PCB (See [Figure 1-16](#) on [page 1-22](#).) and an electronic pressure regulator. The gas serves two functions in the ETD Module:

- As a carrier gas, the nitrogen sweeps the reagent (fluoranthene) from the vial to the ion source where the reagent radical anions are formed.
- As a chemical ionization (CI) vehicle, the nitrogen undergoes collisions with 70 eV electrons from the filament in the ion volume. These 70 eV electrons from the filament knock electrons off of the nitrogen molecules (nitrogen ions are created). The secondary electrons resulting from these collisions have near thermal kinetic energies. These thermal electrons are captured by the fluoranthene to form reagent radical anions that react with the analyte.

Thermo Fisher Scientific strongly recommends a mixture of 25% helium and 75% nitrogen. The helium in this mixture serves as a tracer gas to enable leak checking of gas connections using conventional thermal conductivity-based leak detectors, which are widely used to check leaks in gas chromatography equipment. See [“Gas Supply of the Reagent Ion Source”](#) on [page 1-40](#) for detailed information.

The reagent carrier gas supply in the laboratory is connected to the ETD reagent carrier gas port at the rear side of the ETD Module. See [Figure 1-30](#) on [page 1-40](#).

Reagent Heaters

The reagent heaters (items #9 and #10 in [Figure 1-16](#) on [page 1-22](#), H1 and H2 in [Figure 1-17](#) on [page 1-23](#)) heat the reagent to obtain a sufficient amount of reagent vapor in the carrier gas. The reagent heaters are powered by the Heater Control PCB which, in turn, is controlled by the ETD Control PCB.

The reagent heaters are turned on by selecting the Reagent Ion Source On check box in the Reagent Ion Source dialog box. See [Figure 1-20](#).

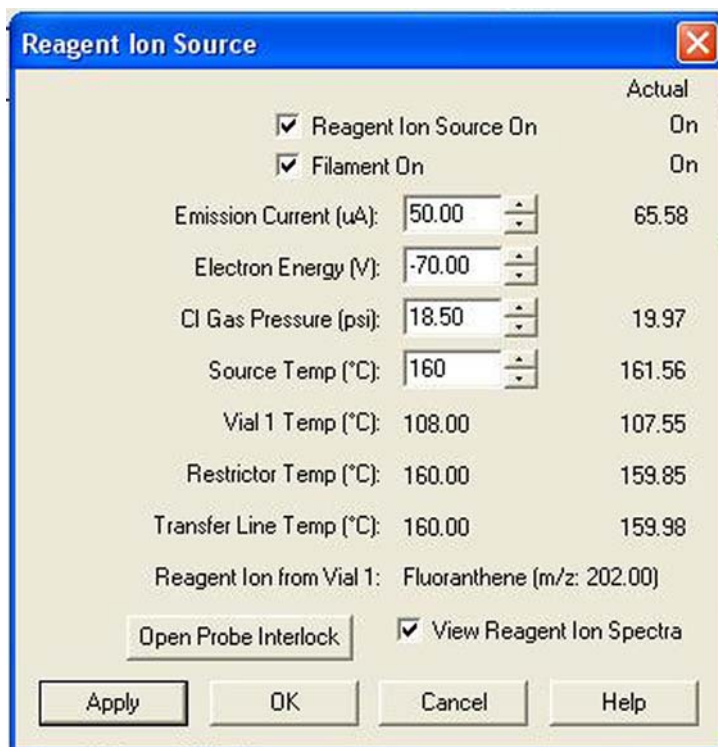


Figure 1-20. Reagent Ion Source dialog box

When you deselect the Reagent Ion Source On check box, the reagent heaters and filament immediately turn off and the reagent ion source goes into Standby mode.

When the reagent ion source is placed in Off mode, cooling nitrogen (high-purity nitrogen) will turn on. This is confirmed by an audible rush (hissing noise) of nitrogen from the reagent ion source area in the back of the ETD Module. This is normal operation. See also [“Turning Off the Reagent Ion Source: What to Expect”](#) on [page 2-14](#).



Warning Burn Hazard. When the reagent ion source is in Off mode, restrictor oven, transfer line, and ion source remain at 160 °C. ▲

The nitrogen cooling gas turns off when the reagent heaters reach 70 °C. If it is necessary to install or replace the reagent vials, follow the procedure in “Replacing the Reagent Vials” on page 3-48.

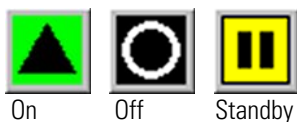
Note The rushing or hissing noise of the nitrogen coming from the back of the ETD Module will stop when the cooling nitrogen turns off. ▲



Warning Burn Hazard. Do not attempt to handle the vials or vial holders when the cooling nitrogen stops. They are still too hot to handle when the cooling nitrogen stops at a vial temperature of 70 °C. Follow the procedures in “Replacing the Reagent Vials” on page 3-48 if it is necessary to install or replace the reagent vials. ▲

When the reagent heaters are in Standby mode, they are immediately turned on by selecting the Reagent Ion Source On check box (Figure 1-20 on page 1-27). When starting from room temperature, it takes up to ten minutes for the reagent heaters and vials to stabilize at 108 °C and reagent to delivered to the ion source.

Note When you switch on a cold reagent ion source, the Tune Plus software warns you that the vial temperature is not sufficient. The filament is automatically switched on after the temperature has stabilized. ▲



When you click the **Standby** button in the Tune Plus window (shown in the left margin), you initiate a Standby process. It delays turning off the reagent heaters and the start of nitrogen cooling for one hour after the system is placed in Standby. This method of placing the system in Standby permits a quick return to operation after a break (lasting one hour or less) rather than waiting up to ten minutes for the heaters to return to temperature and reagent delivery to be fully restored.

In summary:

- The reagent heaters turn off immediately when the Reagent Ion Source On check box is deselected in the Reagent Ion Source dialog box. (See Figure 1-20 on page 1-27.)
- The reagent heaters turn off one hour after the system is placed in Standby by clicking the Standby button in the Tune Plus window.

Reagent Ion Source

The ion source (Figure 1-16 on page 1-22 and inside of the vacuum manifold, see item #14 in Figure 1-16) is where the reagent ions are formed. The ion source contains the filament, the reagent ion volume, and the ion source heater. The filament is the source of electrons that

react with the reagent to form reagent ions. The reagent ion volume is the space where this reaction takes place. See [Figure 1-21](#). The ion source heater is controlled by the ETD Control PCB.

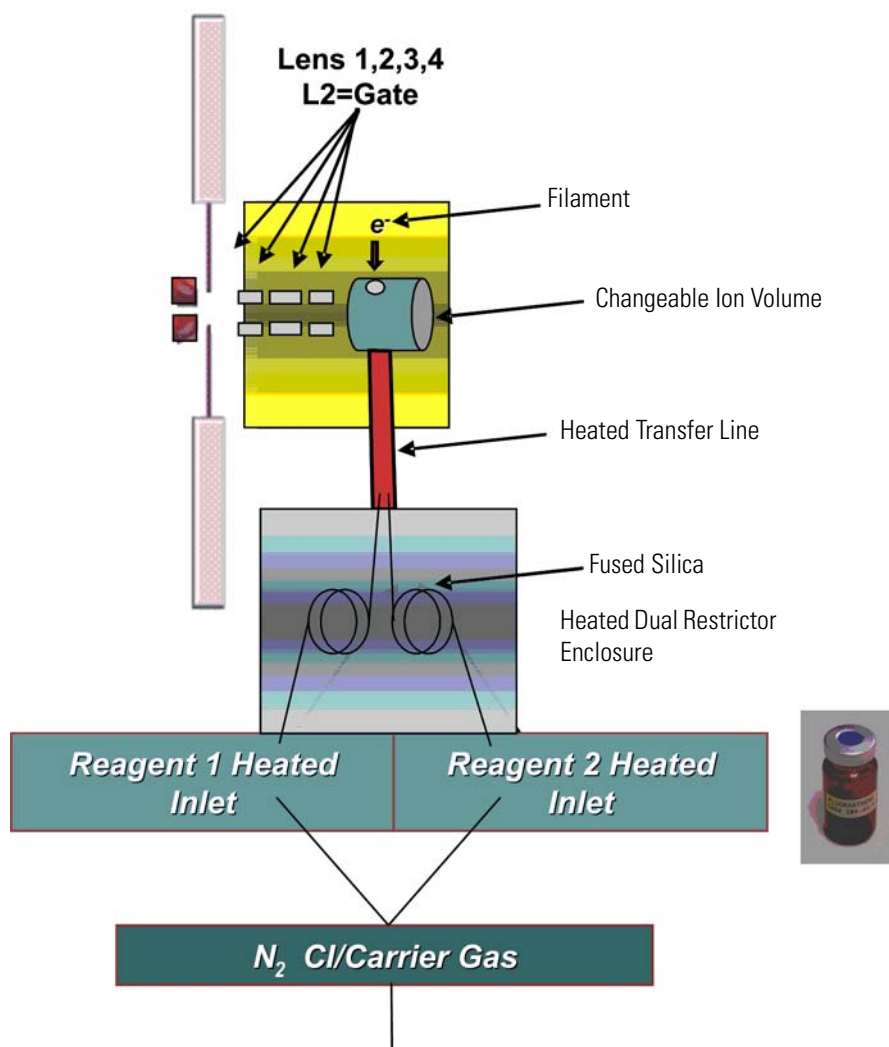


Figure 1-21. Reagent ion source schematics

The reagent ion source contains two reagent vials, CI/carrier gas (nitrogen) handling hardware and flow restrictors, ion volume and filament, ion optics, and heaters for these components. The flow restrictors keep the internal pressure of the reagent vials below atmospheric pressure. This prevents the contents of the reagent vials from being expelled to the laboratory atmosphere.

Vacuum System

Figure 1-22 shows a schematical overview of the Orbitrap analyzer vacuum system.

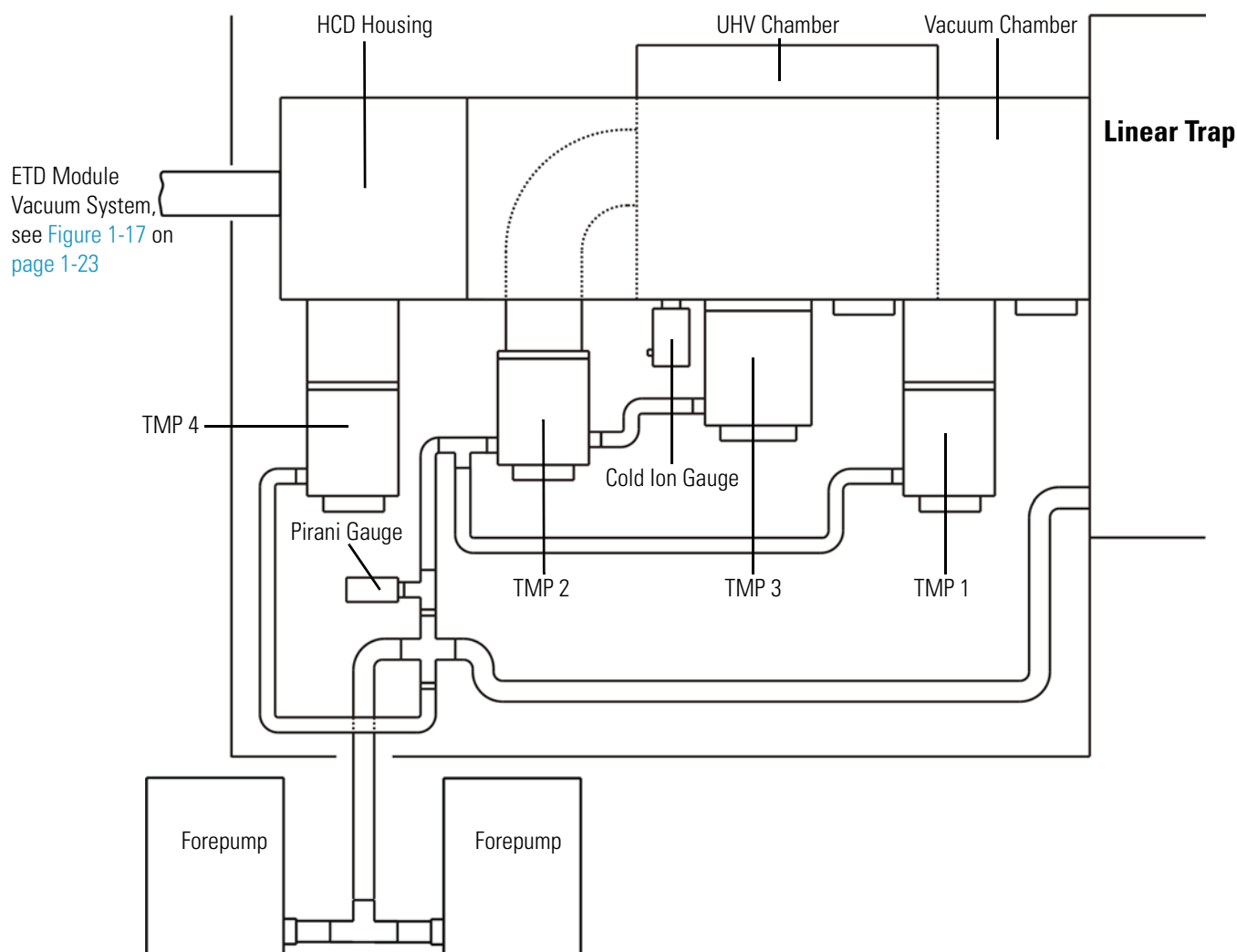


Figure 1-22. Schematic of Orbitrap analyzer vacuum system (CLT compartment and Orbitrap chamber not shown)^a

^aFor an abridged version of the parts list, see [page 4-3](#).

The Orbitrap Velos Pro mass spectrometer has the following vacuum compartments:

- **CLT compartment in the aluminum vacuum chamber** (pumped by the same pump as the linear trap)
- **Vacuum chamber** (pumped by a water-cooled 60 L/s—for N₂—turbomolecular pump HiPace™ 80, *TMP 1*, manufacturer: Pfeiffer)
- **Ultra high vacuum chamber** (UHV chamber, pumped by a water-cooled 60 L/s turbomolecular pump HiPace 80, *TMP 2*, manufacturer: Pfeiffer)

- **Orbitrap chamber** (pumped by a 260 L/s—for N₂—water-cooled turbomolecular pump HiPace 300, *TMP 3*, manufacturer: Pfeiffer)
- **HCD housing** (pumped by a water-cooled 60 L/s turbomolecular pump HiPace 80, *TMP 4*, manufacturer: Pfeiffer)

The forepumps of the linear trap provide the forevacuum for the turbomolecular pumps TMP 1 to TMP 4.

Turbomolecular Pumps

All parts of the system except for the Orbitrap analyzer are mounted in a aluminum vacuum chamber that is evacuated by a 60 L/s turbomolecular pump (TMP 1, see [Figure 1-23](#)). The rotary vane pumps of the linear trap (see [page 1-33](#)) provide the forevacuum for this pump. This chamber is bolted to a stainless steel welded UHV chamber, which accommodates Orbitrap analyzer, lenses, and corresponding electrical connections.

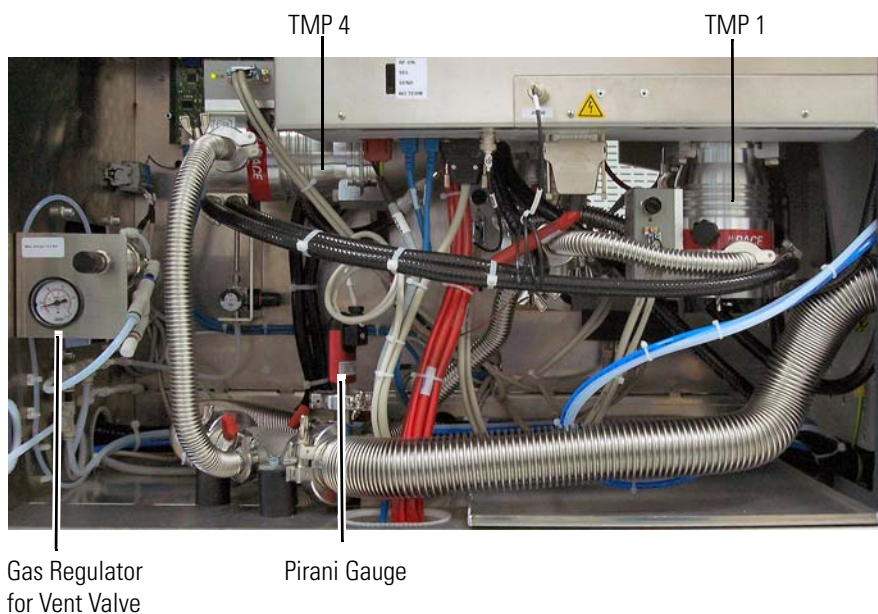


Figure 1-23. Vacuum components on the left instrument side

The UHV chamber is evacuated down to 10⁻⁸ mbar pressure range by a 60 L/s UHV turbomolecular pump (*TMP 2*, see [Figure 1-24](#) on [page 1-32](#)). The Orbitrap analyzer itself is separated from the UHV chamber by differential apertures and is evacuated down to 10⁻¹⁰ mbar by a 260 L/s turbomolecular pump (*TMP 3*, see [Figure 1-24](#)). The HCD housing is evacuated by a 60 L/s UHV turbomolecular pump (*TMP 4*, see [Figure 1-23](#)) that is mounted to its bottom via an elbow. This dedicated pump for the HCD cell protects the low pressure cell of the linear trap from gas overload.

In the Orbitrap Velos Pro ETD mass spectrometer, a tube that contains the transfer multipole (flatapole) connects the HCD housing to the ETD Module.

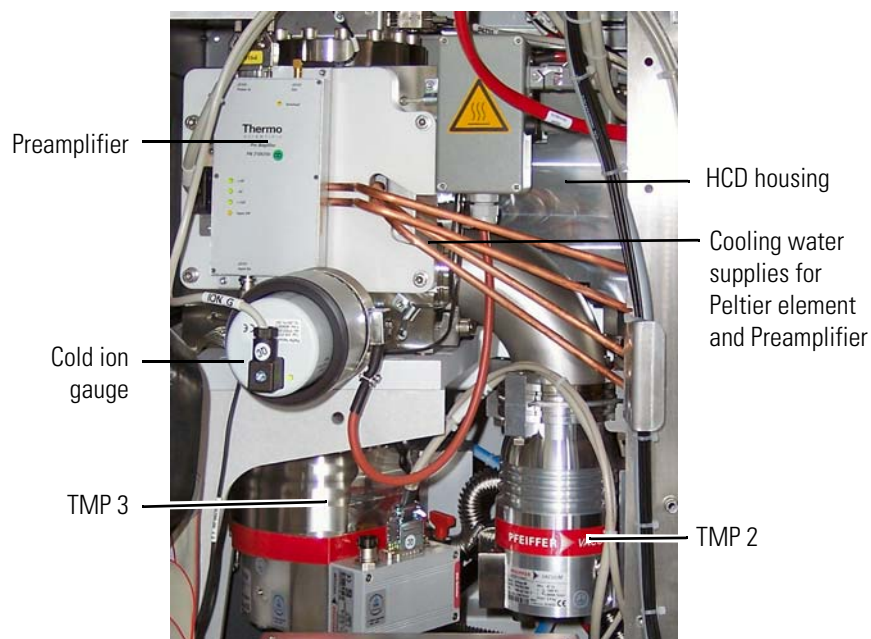


Figure 1-24. Vacuum components on the right instrument side

All turbomolecular pumps are equipped with TC 110 control units (manufacturer: Pfeiffer). A 24 V switch mode power supply provides the electric power for all four turbomolecular pumps of the system.

Linear Trap Turbomolecular Pump

A separate turbomolecular pump provides the high vacuum for the linear ion trap. It is mounted to the bottom of the vacuum manifold of the linear ion trap. For more information, refer to the *LTQ Series Hardware Manual*.

ETD Module Turbomolecular Pump

In the Orbitrap Velos Pro ETD mass spectrometer, a separate turbomolecular pump (Edwards EXT75DX) provides the high vacuum for the ETD reagent ion source. See [Figure 1-16](#) on [page 1-22](#). It is backed up by a dedicated rotary vane pump at the bottom of the ETD Module. See [Figure 1-26](#) on [page 1-34](#). This air-cooled turbomolecular pump contains no user-serviceable parts.

Forepumps of the Linear Trap

The rotary vane pumps from the linear trap serve as forepumps for the three smaller turbomolecular pumps (TMP 1, TMP 2, and TMP 4). The exhaust hose from the forepumps is led to the back of the instrument and connects them to the exhaust system in the laboratory. The forepumps are located on a small cart in the forepumps cabinet below the linear trap. See [Figure 1-25](#).

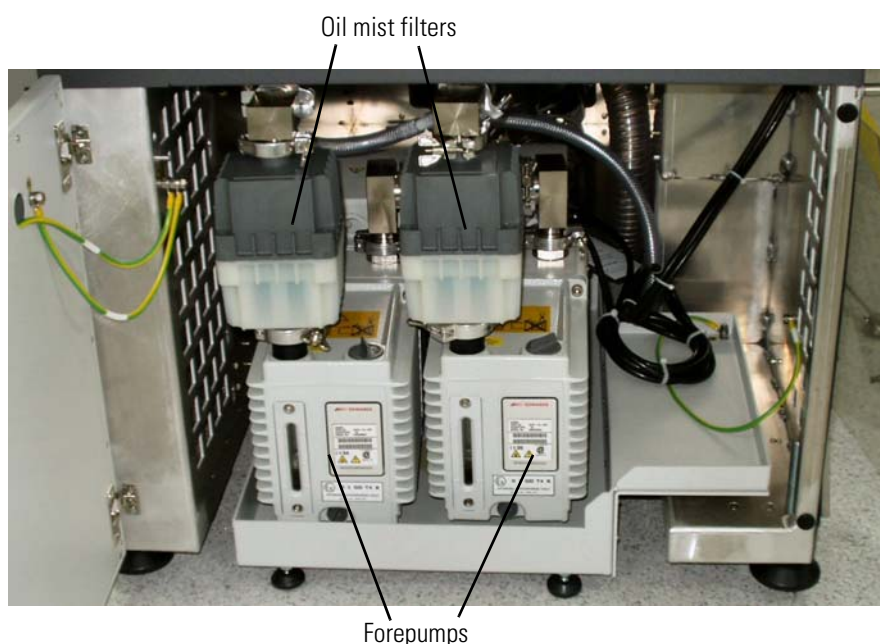


Figure 1-25. Forepumps cabinet

To minimize the ingress of pump oil into the exhaust system, the outlets of the forepumps are fitted to oil mist filters. See [page 3-11](#) on instructions about returning the collected oil to the forepumps.

The forepumps of the linear trap are powered by the power panel of the linear ion trap.



Warning Health Hazard. Hazardous materials may be present in the effluent of the forepumps. The connection to an adequate exhaust system is mandatory! ▲

Leave the switches of the forepumps always in the On position to provide the control from the vacuum control panel. Before starting the pumps, however, make sure that:

- The forevacuum pumps are filled with oil,
- They are connected to the power supply, and
- The gas ballast is shut.

For a detailed description of the forepumps, refer to the handbook of the manufacturer.

Forepump of the ETD Module

In the Orbitrap Velos Pro ETD mass spectrometer, a rotary vane pump (Edwards RV 3) provides the forevacuum for the ETD turbomolecular pump. (See “[ETD Module Turbomolecular Pump](#)” on [page 1-32](#).) It is located in a cabinet at the bottom of the ETD Module. The ETD forepump is equipped with an oil mist filter and stands on a drip pan. See [Figure 1-26](#) on [page 1-34](#).

An exhaust hose connects the forepump to the exhaust system in the laboratory. A forevacuum tube connects the ETD forepump to the ETD TMP. The forepump electrical cord is plugged into the Forepump receptacle on the ETD Module power panel. See [Figure 1-19](#) on [page 1-24](#).

For maintenance instructions for the ETD forepump, see “[Maintenance of the ETD Forepump](#)” on [page 3-5](#) and the manual that came with the forepump.

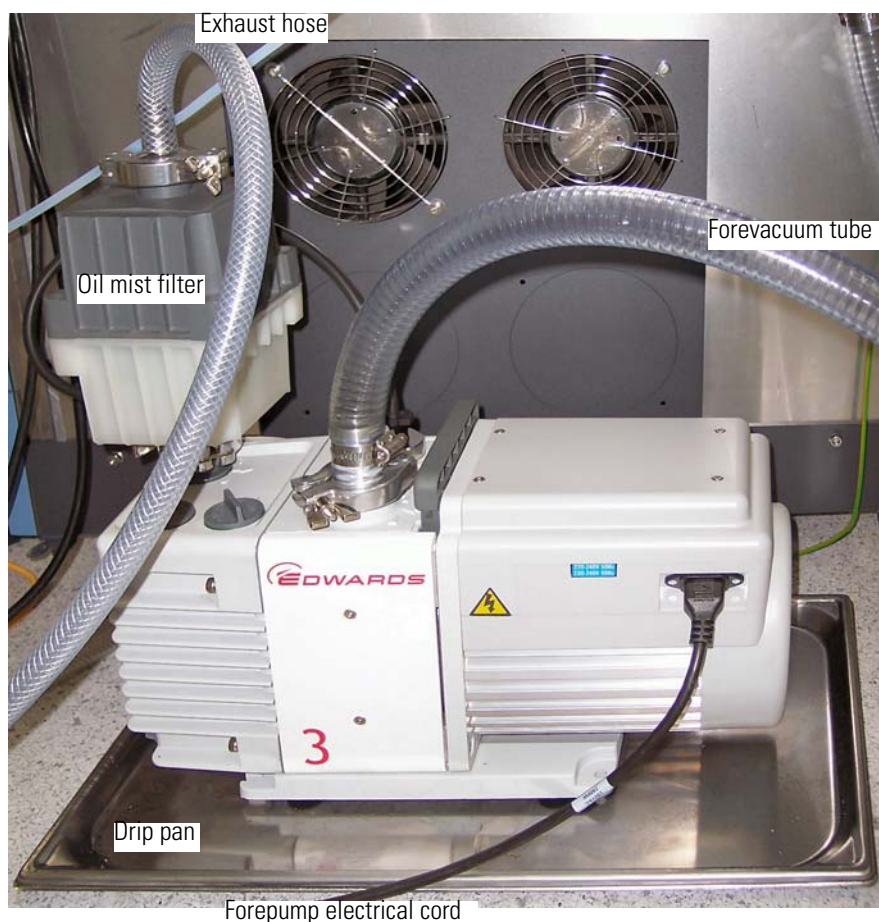


Figure 1-26. Forepump for ETD Module

Vacuum System Controls

The power distribution board controls all turbomolecular pumps via voltage levels. See “[Power Distribution Board](#)” on [page 1-53](#). An interface for RS485 data via the instrument control board connects the turbomolecular pumps with the linear ion trap. (See “[Instrument Control Board](#)” on [page 1-52](#).) The turbomolecular pump of the linear ion trap and the ETD turbomolecular pump have individual controllers.

Vacuum Gauges

Several vacuum gauges monitor the vacuum within the instrument:

- The forevacuum of the Orbitrap Velos Pro mass spectrometer is monitored by an Active Pirani gauge (TPR 280, manufacturer: Pfeiffer) connected to the forevacuum line. See [Figure 1-23](#) on [page 1-31](#).
- The high vacuum of the Orbitrap Velos Pro mass spectrometer is monitored by a Cold Ion Gauge (IKR 270, manufacturer: Pfeiffer) connected to the UHV chamber. See [Figure 1-24](#) on [page 1-32](#). Because the gauge would be contaminated at higher pressures, it is turned on only when the forevacuum has fallen below a safety threshold ($<10^{-2}$ mbar).
- The linear ion trap vacuum is monitored by a Convectron™ gauge and an ion gauge. Refer to the *LTQ Series Hardware Manual* for more information.
- In the Orbitrap Velos Pro ETD mass spectrometer, two dedicated vacuum gauges monitor the vacuum in the ETD Module. A Convectron gauge (see [Figure 1-16](#) on [page 1-22](#) and [Figure 1-17](#) on [page 1-23](#)) monitors the pressure in the ETD forevacuum line and an ion gauge (see [Figure 1-16](#)) monitors the pressure in the reagent ion source. [Table 1-3](#) shows typical pressure readings in the ETD Module.

Table 1-3. Typical pressure readings in the ETD Module

Conditions	Convection Gauge Reading	Ion Gauge Reading
CI gas pressure set to 20 psi	0.1–0.01 Torr	$20\text{--}35 \times 10^{-5}$ Torr

The vacuum gauges of the Orbitrap Velos Pro mass spectrometer are connected to the power distribution board that directly responds to the pressure values. (See “[Power Distribution Board](#)” on [page 1-53](#).) The analog values are digitized by the instrument control board. (See “[Instrument Control Board](#)” on [page 1-52](#).) They are then sent as readout values to the data system.

Switching on the Vacuum System

When the vacuum system is switched on, the following occurs:

1. After the Vacuum Pumps switch is switched On, the pumps of the linear ion trap and the Orbitrap Velos Pro mass spectrometer are run up. The Pirani gauge (see above) controls the Orbitrap Velos Pro MS low vacuum pressure as well as the pressure at the forevacuum pumps. Within a short time, a significant pressure decrease must be observed. The goodness of the vacuum can be estimated by means of the rotation speed of the turbomolecular pumps (for example, 80% after 15 minutes).
2. If the working pressure is not reached after the preset time, the complete system is switched off. At the status LED panel of the power distribution board, an error message (Vacuum Failure) is put out (see below).
3. The Cold Ion Gauge is only switched on after the low vacuum is reached. It is then used to monitor the vacuum in the Orbitrap analyzer region.

Vacuum Failure

In case the pressure in the Orbitrap Velos Pro mass spectrometer or the linear ion trap exceeds a safety threshold, the complete system including linear ion trap, electronics, and pumps is switched off. However, the power distribution is kept under current and puts out an error message at the LED panel. (See [“Power Distribution Board”](#) on [page 1-53](#).) It can be reset by switching the main power switch off and on. (See [“Main Power Switch”](#) on [page 1-9](#).)

Upon venting, the vent valves of the turbomolecular pumps on the Orbitrap analyzer stay closed. Only the vent valve of the linear ion trap is used. (See [“Vent Valve of the Linear Ion Trap”](#) on [page 1-39](#).)

Vacuum System Heating during a System Bakeout

After the system has been open to the atmosphere (for example, for maintenance work), the vacuum deteriorates due to contaminations of the inner parts of the vacuum system caused by moisture or a power outage. These contaminations must be removed by heating the vacuum system: a system bakeout. See [“Baking Out the System”](#) on [page 3-4](#).

Gas Supply

This section describes the gas supplies for the mass analyzers of the Orbitrap Velos Pro mass spectrometer and the reagent ion source of the Orbitrap Velos Pro ETD mass spectrometer.

Gas Supply for the Mass Analyzers

Figure 1-27 shows a schematical view of the gas supply for the instrument. The gas supply of the ETD system is highlighted in gray.

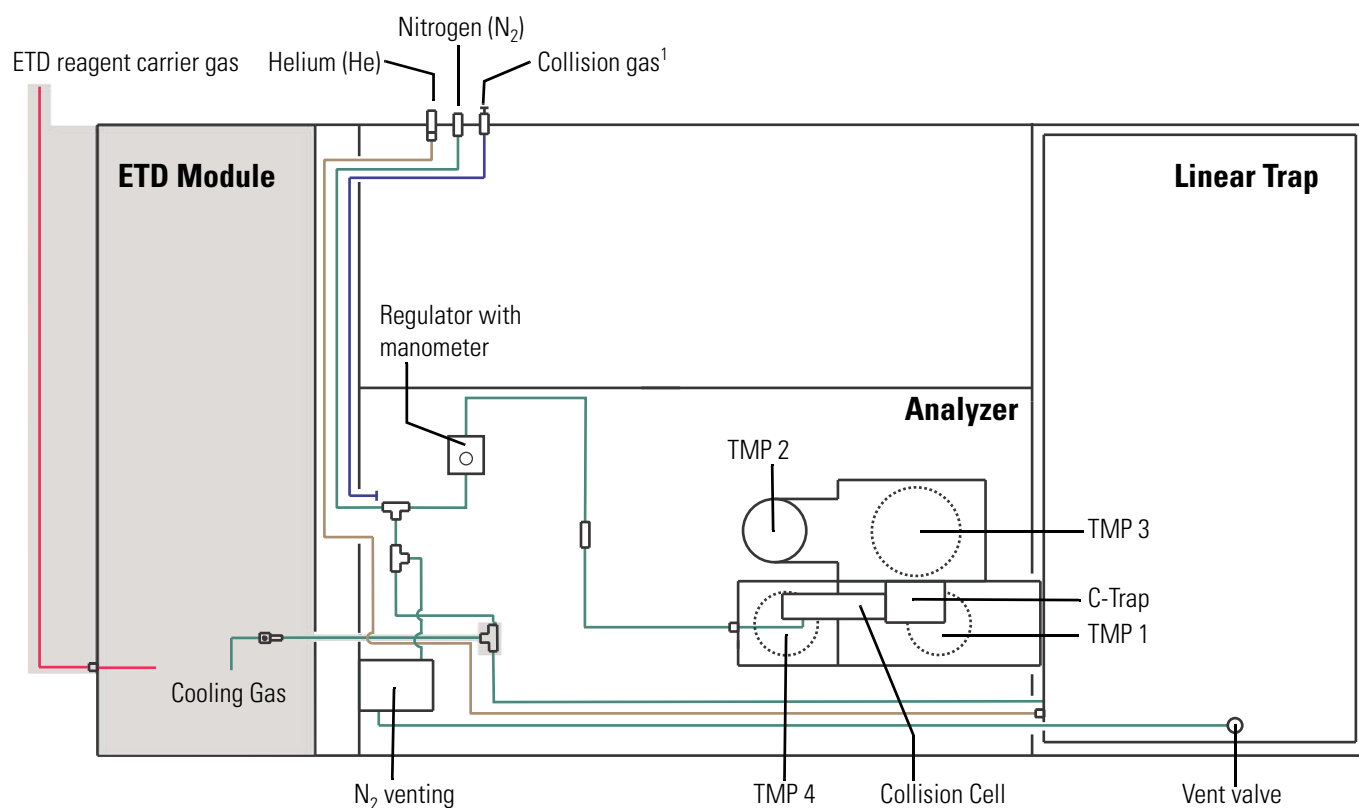


Figure 1-27. Schematic of gas supply for Orbitrap Velos Pro ETD MS^a

^aFor parts lists of the gas supply, see [page 4-4](#) and [page 4-5](#).

Gas Inlet Ports of the Instrument

On its right side (See [Figure 1-9](#) on [page 1-10](#).), the instrument provides three gas inlet ports for the gas supply of the mass analyzers:

- Nitrogen: The linear trap requires high-purity (99%) nitrogen for the API sheath gas and auxiliary/sweep gas. The required gas pressure is 690 ± 140 kPa (6.9 ± 1.4 bar, 100 ± 20 psi).

¹ The port named Collision Gas is not used in the Orbitrap Velos Pro MS.

In the Orbitrap Velos Pro ETD mass spectrometer, the ETD system uses the high-purity nitrogen for cooling the reagent vials when the reagent ion source is turned off.

- Helium: The linear trap requires ultra-high purity (99.999%) helium for the collision gas. The required gas pressure is 275 ± 70 kPa (2.75 ± 0.7 bar, 40 ± 10 psi).



Warning Danger of Asphyxiation. Accumulation of nitrogen gas could displace sufficient oxygen to suffocate personnel in the laboratory. Ensure that the laboratory is well ventilated. ▲

❖ **To connect the nitrogen source to the Orbitrap Velos Pro mass spectrometer**

1. Connect an appropriate length of Teflon™ tubing to the nitrogen source in the laboratory. The Installation Kit contains 6 m (20 ft) of suitable Teflon tubing (OD 6 mm, P/N 0690280). The connection for the Teflon hose to the nitrogen gas supply is not provided in the kit; you have to supply this part.
2. Connect the opposite end of the Teflon tubing to the press-in fitting labeled Nitrogen, which is located at the right side of the instrument. See [Figure 1-9](#) on [page 1-10](#). To connect the tubing, align the Teflon tubing with the opening in the fitting and firmly push the tubing into the fitting until the tubing is secure.

❖ **To connect the helium source to the Orbitrap Velos Pro mass spectrometer**

1. Connect an appropriate length of 1/8-in. ID copper or stainless steel tubing with a brass Swagelok-type 1/8-in. nut (P/N 00101-15500) and a 2-piece brass 1/8-in. ID ferrule [P/N 00101-08500 (front), P/N 00101-2500 (back)] to the Helium gas inlet. See [Figure 1-28](#) for the proper orientation of the fitting and ferrule.
2. Connect the opposite end of the tubing to the helium gas source, using an appropriate fitting.

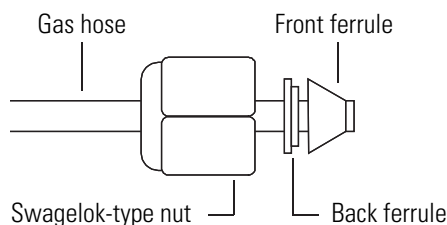


Figure 1-28. Proper orientation of the Swagelok-type nut and two-piece ferrule

Gas Distribution Within the Instrument

Helium gas is led from the helium port through a stainless steel capillary to the right rear side of the linear trap. See [Figure 1-27](#) on [page 1-37](#). High purity nitrogen gas is led from the nitrogen port via Teflon tubing to the right side of the Orbitrap Velos Pro mass spectrometer. Here, two T-pieces divide the nitrogen gas flow into three parts.

The first part of the high purity nitrogen gas flow is directed through Teflon tubing via a pressure regulator to the vent valve of the linear trap. (See [Figure 1-29](#) on [page 1-39](#).) The second part of the nitrogen flow is directed through Teflon tubing to the API source. The third part of the nitrogen flow enters a gas regulator with manometer, which keeps the gas pressure to the C-Trap and HCD collision cell constant. (See [Figure 1-29](#), background.) From the regulator, the collision gas is led through *red* PEEKSil™ tubing (100 mm ID silica capillary in 1/16 inch PEEK tubing) to the collision octapole next to the curved linear trap (flow rate: ~0.5 mL/min). The nitrogen gas leaking from the HCD collision cell (3–5 mbar) is used for ion trapping and cooling in the C-Trap.

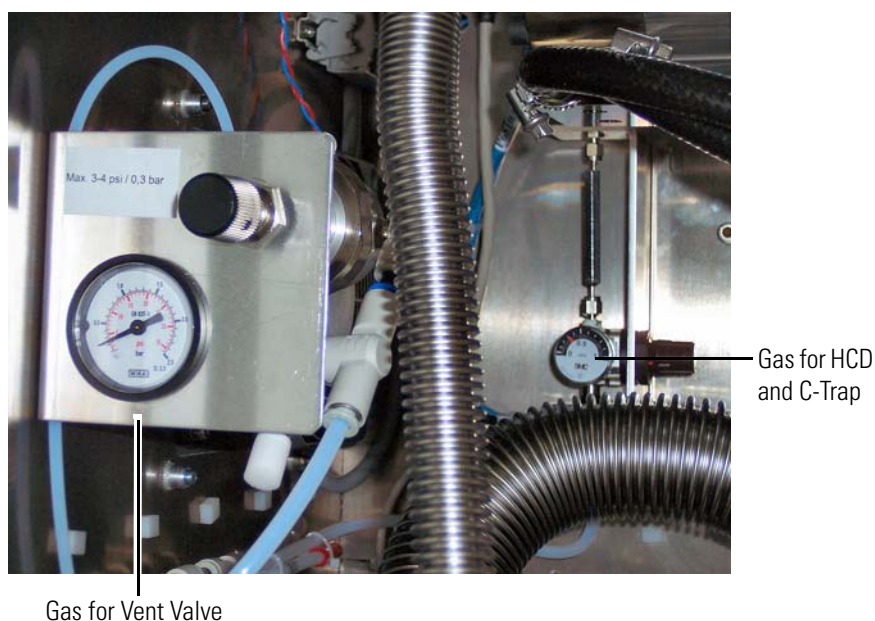


Figure 1-29. Gas regulators

In the Orbitrap Velos Pro ETD mass spectrometer, nitrogen is also directed through Teflon tubing to the ETD Module to be used for cooling the reagent vials when the reagent ion source is turned off.

Vent Valve of the Linear Ion Trap

If the system and pumps are switched off, the system is vented. The vent valve is controlled by the linear ion trap. The *LTQ Series Hardware Manual* contains further information about the vent valve.

The instrument is vented with high purity nitrogen from the same tubing that supplies the Velos Pro MS sheath gas. See [Figure 1-27](#) on [page 1-37](#). The vent valve of the Velos Pro mass spectrometer is supplied via a pressure regulator that is set to a venting pressure of 3–4 psi. The pressure regulator is located at the left side of the Orbitrap Velos Pro mass spectrometer. (See [Figure 1-29](#), front.)

Gas Supply of the Reagent Ion Source

In addition to high purity nitrogen for cooling, the reagent ion source of the Orbitrap Velos Pro ETD mass spectrometer uses a mixture of 25% helium and 75% nitrogen gas as carrier gas and chemical ionization (CI) vehicle. This gas mixture must be ultra high-purity (minimum purity 99.999%) with less than 3.0 ppm each of water, oxygen, and total hydrocarbons. The required gas pressure is 690 ± 140 kPa (6.9 ± 1.4 bar, 100 ± 20 psi). The ETD carrier gas supply of the laboratory is connected via metal tubing to the inlet port at the rear side of the instrument. See [Figure 1-30](#).



Figure 1-30. ETD reagent carrier gas port at the ETD Module

The helium in this mixture serves as a tracer gas to enable leak checking of gas connections using conventional thermal conductivity-based leak detectors, which are widely used to check leaks in gas chromatography equipment.

Note If the helium/nitrogen mixture is not available, then use a nitrogen supply that is ultra high-purity (99.999%) with less than 3.0 ppm each of water, oxygen, and total hydrocarbons. ▲

Triple Gas Filter

A triple (oxygen/water/hydrogen) gas filter is installed between the regulator on the reagent carrier gas source and the ETD module to ensure that the reagent carrier gas (either nitrogen or helium/nitrogen) is better than 99.999% pure with much less than 1 ppm of oxygen, water, and hydrocarbons.

Refer to the filter manufacturer's instructions for information about how to monitor the color changes in the filters that indicated when the filters need to be replaced, as well as information about where to order new filters. If there are no leaks in the reagent carrier gas plumbing, you can expect the filters to last a year or more. Thermo Fisher Scientific strongly recommends that a Thermo Fisher Scientific field service engineer replace the gas filters.

Cooling Water Circuit

Figure 1-31 on page 1-42 shows a schematical view of the cooling water circuit in the Orbitrap Velos Pro mass spectrometer. For a parts list of the cooling water circuit, see page 4-4. Cooling water at a temperature of 20 °C enters and leaves the instrument at the bottom of the right side. See Figure 1-9 on page 1-10. First, the fresh water passes through the turbomolecular pumps in the order TMP 3 → TMP 1 → TMP 4 → TMP 2. Then it passes through the heating element (Peltier element) that maintains (± 0.5 °C) the preset temperature of the analyzer. After that, the cooling water passes through the preamplifier cooling unit. Before it leaves the instrument, the water passes through the other Peltier element at the back of the central electrode power supply board.

A flow control sensor is connected to the power distribution board and allows displaying the current flow rate of the cooling water in the software. An inline filter, which is installed upstream, protects the sensor. It must be replaced annually, see page 3-58 for instructions.

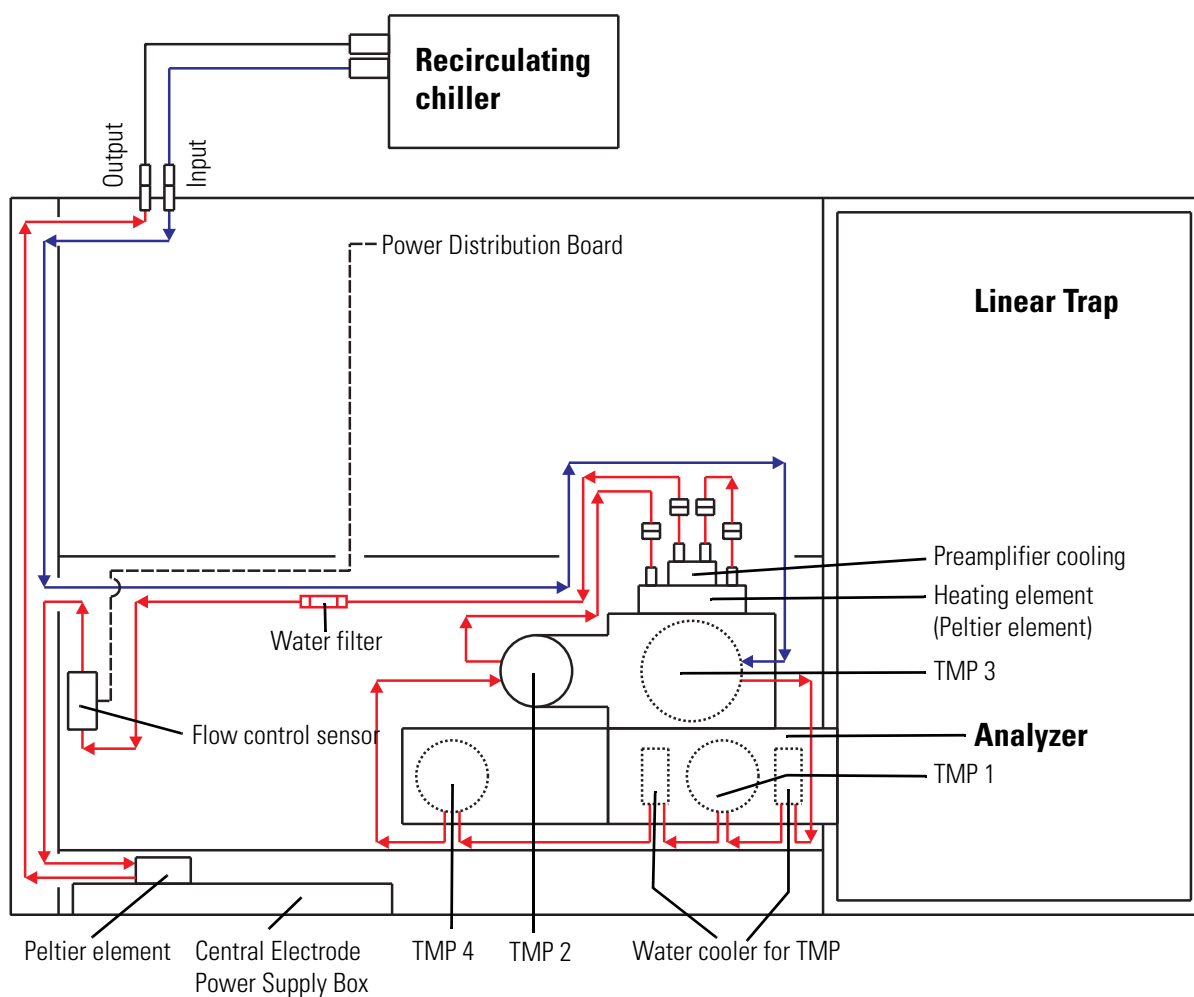


Figure 1-31. Schematic of cooling water circuit

Recirculating Chiller

A recirculating chiller (Thermo Scientific NESLAB ThermoFlex™ 900) is shipped with the instrument, making the mass spectrometer independent from any cooling water supply. A wall receptacle provides the electric power for the chiller. Two water hoses (black), internal diameter 9 mm, wall thickness 3 mm, length approx. 3 m (~10 ft) are shipped with the instrument.

For instruction about performing maintenance for the chiller, see “Maintenance of the Cooling Circuit” on page 3-58. See also the manufacturer’s manual for the chiller.

Properties of Cooling Water

The water temperature is not critical, but should be in the range of 20 to 25 °C (68 to 77 °F). Lower temperatures could lead to a condensation of atmospheric water vapor. It is recommended to use distilled water rather than de-ionized water due to lower concentration of bacteria and residual organic matter.

The water should be free of suspended matter to avoid clogging of the cooling circuit. In special cases, an in-line filter is recommended to guarantee consistent water quality.

The cooling water should meet the following requirements:

Hardness:	<0.05 ppm
Resistivity:	1–3 MΩ/cm
Total dissolved solids:	<10 ppm
pH:	7–8



Warning Burn Hazard. If the water circuit fails, all parts of the water distribution unit may be considerably heated up. Do not touch the parts! Before disconnecting the cooling water hoses, make sure the cooling water has cooled down! ▲

Printed Circuit Boards

The Orbitrap Velos Pro mass spectrometer is controlled by a PC running the Xcalibur™ software suite. The software controls all aspects of the instrument. The main software elements are the communication with the linear ion trap, the control of ion detection, and the control of the Orbitrap analyzer.

The following pages contain a short overview of the electronic boards in the MS portion of the Orbitrap Velos Pro mass spectrometer. For each board, its respective location and function are given. If applicable, the diagnostic LEDs on the board are described. For a description of the printed circuit boards in the ETD Module, see “[ETD Module](#)” on [page 1-21](#).

The electronics of the Orbitrap Velos Pro mass spectrometer contains complicated and numerous circuits. Therefore, only qualified and skilled electronics engineers should perform servicing.

A Thermo Fisher Scientific field service engineer should be called if servicing is required. It is further recommended to use Thermo Fisher Scientific spare parts only. When replacing fuses, only use the correct type. Before calling a service engineer, please try to localize the defect via errors indicated in the software or diagnostics. A precise description of the defect will ease the repair and reduce the costs.



Warning Electrical Shock Hazard. Parts of the printed circuit boards are at high voltage. Shut down the instrument and disconnect it from line power before performing service. Opening the electronics cabinet is only allowed for maintenance purposes by qualified personnel. ▲

Note Many of the electronic components can be tested by the Orbitrap Velos Pro MS diagnostics, which is accessible from the Tune Plus window. ▲

Linear Ion Trap Electronics

The linear ion trap is connected to the Orbitrap Velos Pro MS main power switch. The linear ion trap has a sheet metal back cover.

[Figure 1-32](#) shows the electronic connections at the rear side of the linear trap.



Figure 1-32. Electronic connections to linear trap

The linear ion trap electronics has two connections with the Orbitrap Velos Pro MS electronics:

- Data communication with the internal computer of the Orbitrap Velos Pro mass spectrometer. See [“Electronic Boards on the Right Side of the Instrument”](#) on [page 1-46](#).
- Signal communication (SPI bus) with supply information for the instrument control board. See [“Instrument Control Board”](#) on [page 1-52](#).

For further information about the linear ion trap electronics, refer to the *LTQ Series Hardware Manual*.

Electronic Boards on the Right Side of the Instrument

Figure 1-33 shows the parts of the instrument when the right side panel is opened. A transparent cover protects the lower part.

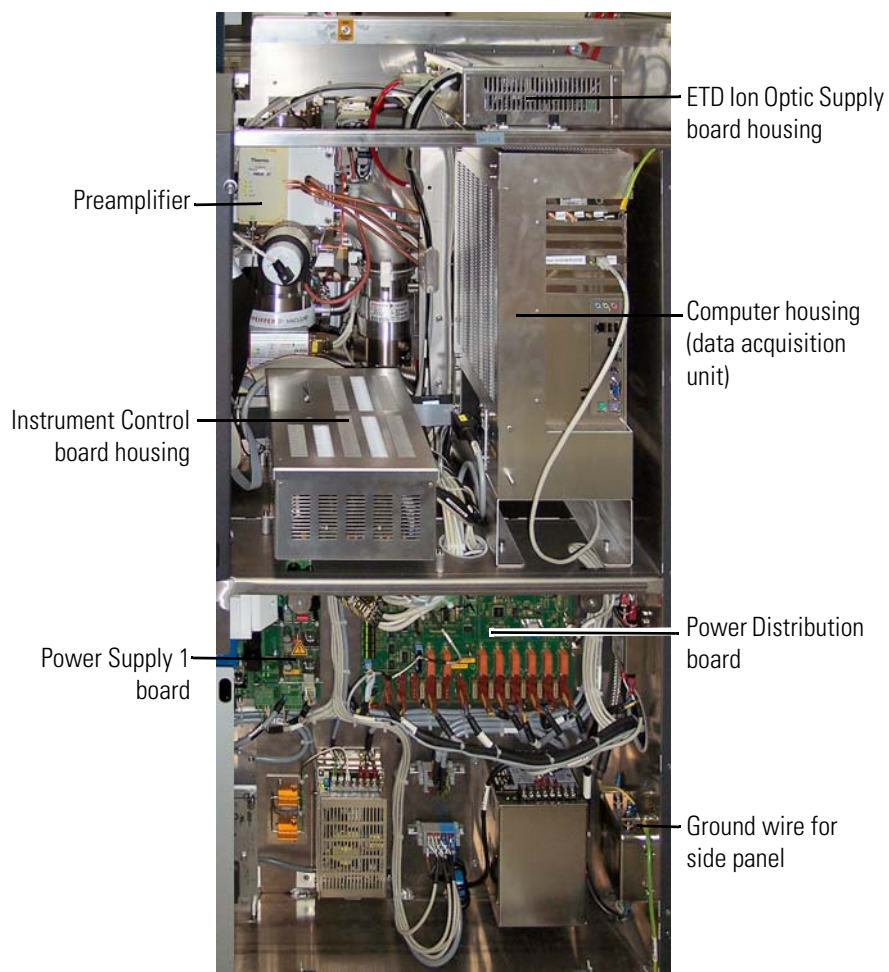


Figure 1-33. Electronic boards on the right side of the instrument



The side panel is connected to the instrument frame by two green/yellow ground wires. See bottom of Figure 1-33. The connectors on the panel are labeled with green-yellow PE (for **P**rotective **E**arth) signs. See photo left. Do not forget to reconnect them before closing the panel!

ETD Ion Optic Supply Board

The ETD Ion Optic Supply board is mounted on top of the data acquisition unit. See Figure 1-34. It supplies the voltages for the HCD collision cell. In the Orbitrap Velos Pro ETD mass spectrometer, this board also supplies the RF voltage and the DC voltages for the ETD Module: an RF voltage with DC offset, three DC voltages with ± 250 V, and a DC voltage with ± 12 V.

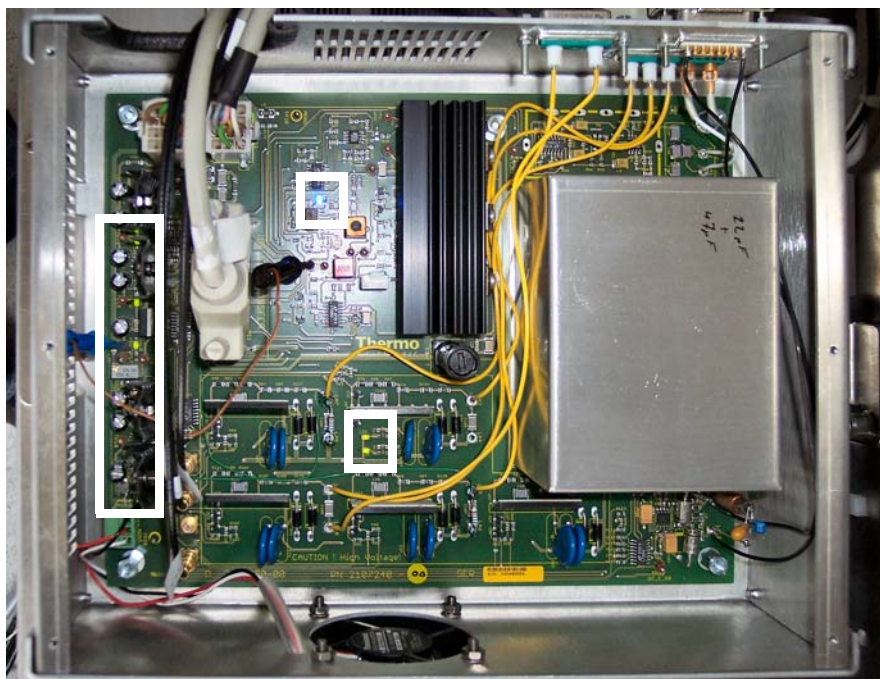


Figure 1-34. ETD Ion Optic Supply board

The diagnostic LEDs on the ETD ion optic supply board are listed in [Table 1-4](#) on [page 1-47](#). The positions of the diagnostic LEDs on the board are indicated by white rectangles in [Figure 1-35](#).

Table 1-4. Diagnostic LEDs on the ETD Ion Optic Supply board

No.	Name	Color	Description	Normal Operating Condition
LD1	+275 V	Green	+275 V input voltage present	On
LD2	-275 V	Green	-275 V input voltage present	On
LD3	RF Supply	Green	RF input voltage (22 V) present	On
LD4	+24 V	Green	+24 V input voltage present	On
LD5	+15 V	Green	+15 V input voltage present	On
LD6	-15 V	Green	-15 V input voltage present	On
LD7	RF1_ON	Blue	RF-generator switched on	On/Off, depending on active application

Preamplifier

The preamplifier is located in a housing next to the Cold Ion Gauge. See [Figure 1-35](#). It is water cooled to protect it during a system bakeout.

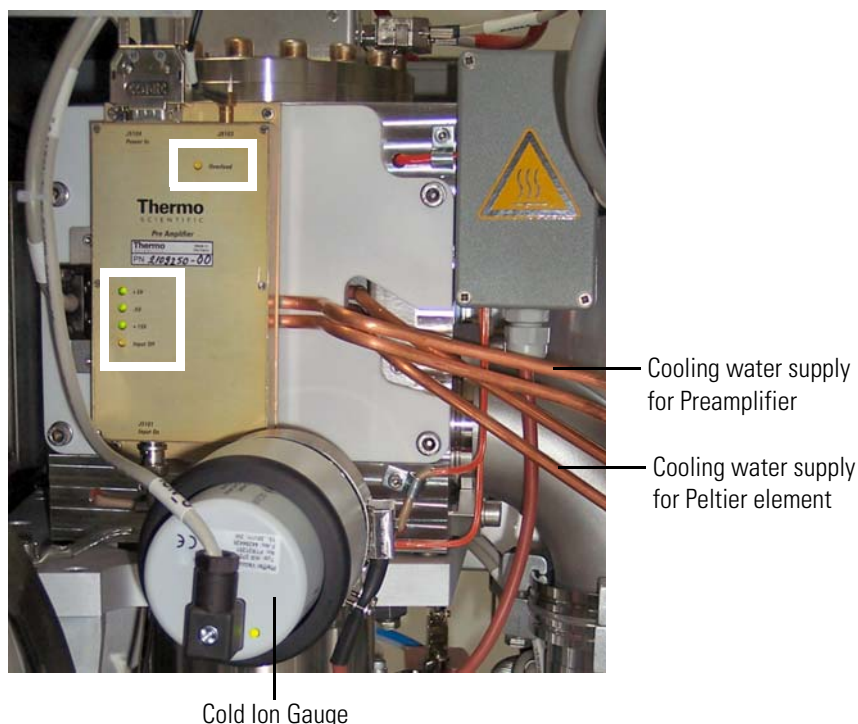


Figure 1-35. Preamplifier board

This board is a broadband preamplifier with differential high-impedance inputs. It serves as a detection amplifier and impedance converter for the image current created by the oscillating ions. The output current is transferred to the data acquisition board. It has an amplification factor of about 60 dB and covers the frequency range from 15 kHz to 10 MHz.

The diagnostic LEDs on the preamplifier are listed in [Table 1-5](#) on [page 1-48](#). The positions of the diagnostic LEDs on the board are indicated by white rectangles in [Figure 1-35](#).

Table 1-5. Diagnostic LEDs on the Preamplifier board

No.	Name	Color	Description	Normal Operating Condition
LD1	Overload	Yellow	RF output is overloaded	Off
LD2	+5 V	Green	+5 V input voltage present	On
LD3	+15 V	Green	+15 V input voltage present	On
LD4	-5 V	Green	-5 V input voltage present	On
LD5	Input off	Yellow	RF inputs are shortened (protection)	On, off during Detect

Internal Computer

Figure 1-36 shows the components of the data acquisition unit. The unit is mounted in a housing located at the right side of the instrument.

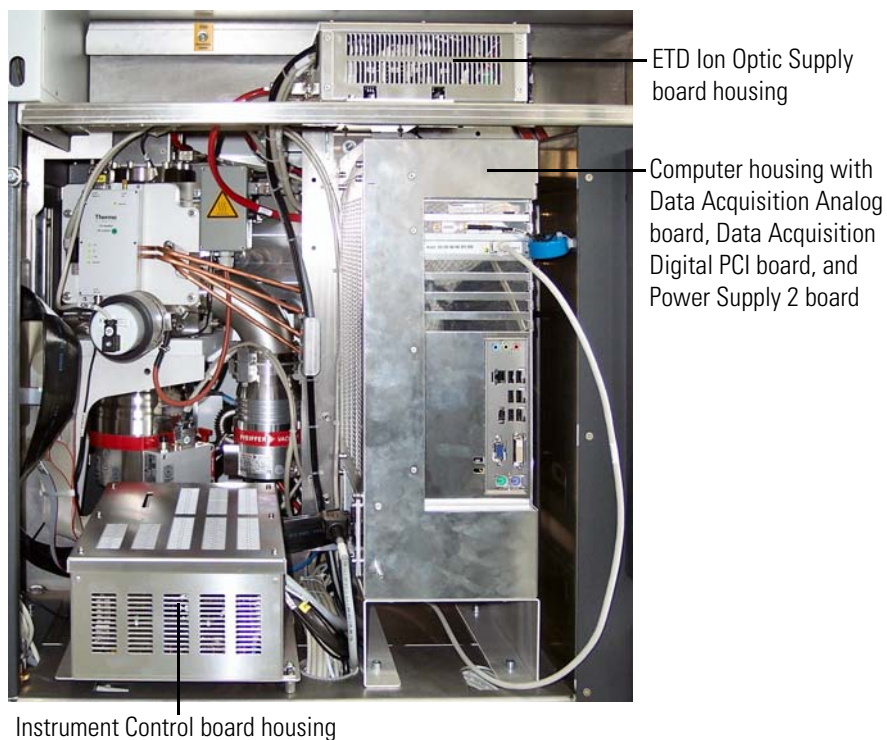


Figure 1-36. Data Acquisition unit

The internal computer contains a computer mainboard with an ATX power supply. The data acquisition digital PCI board is directly plugged into the mainboard. The data acquisition analog board is mounted on top of the computer mainboard.

Data Acquisition Digital PCI Board

Figure 1-37 shows the data acquisition digital PCI board. It is an add-on board to the internal computer. (See Figure 1-36 on page 1-49.)

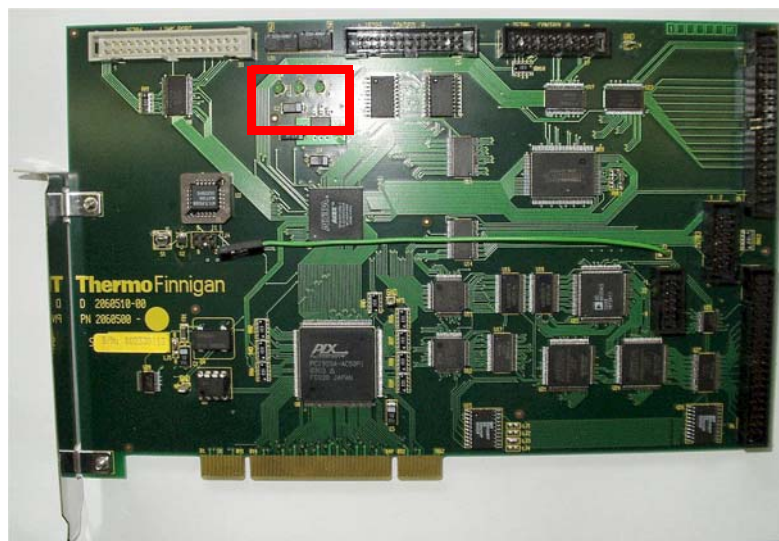


Figure 1-37. Data Acquisition Digital PCI board

This board is used to convert detected ion signals to digital form and to interface to the computer mainboard. The board has two 16 bit parallel connections to the DAC and the ADC on the data acquisition analog board, which are used for controlling and reading-back signals. A high-speed link port channel is also on the board that is used to communicate with the electronics in the linear ion trap.

Precision timing is derived from the data acquisition analog board and events with lower requirements use the timer in the internal computer. This timer is used to check at regular intervals whether the foreground process works as expected.

Communication takes place not only between the linear ion trap and the internal computer of the Orbitrap Velos Pro system, but also between the linear ion trap and the data system computer. For further information about the data system, refer to the *LTQ Series Hardware Manual*.

The diagnostic LEDs listed in Table 1-6 show the status of the board. The position of the LEDs on the board is indicated by a red rectangle in Figure 1-37.

Table 1-6. Diagnostic LEDs of the Data Acquisition Digital PCI board

Name	Color	Description	Normal Operating Condition
+5 V	Green	+5 V voltage present	On
+3.3 V	Green	+3.3 V voltage present	On
+2.5 V	Green	+2.5 V voltage present	On

Data Acquisition Analog Board

Figure 1-38 shows the data acquisition analog board. This board is an add-on board to the mainboard of the internal computer. See Figure 1-36 on page 1-49. It is used to convert analog to digital signals for Orbitrap analyzer experiments, especially for detecting the ions. The board contains an ADC for the detection of the transient signal, with a frequency range from 10 kHz to 10 MHz. Three anti-aliasing filters for the low, middle and high mass range are automatically selected by the software.

The data acquisition board provides precision timing to control the acquisition. Events with lower timing requirements on accuracy are controlled by the linear ion trap.

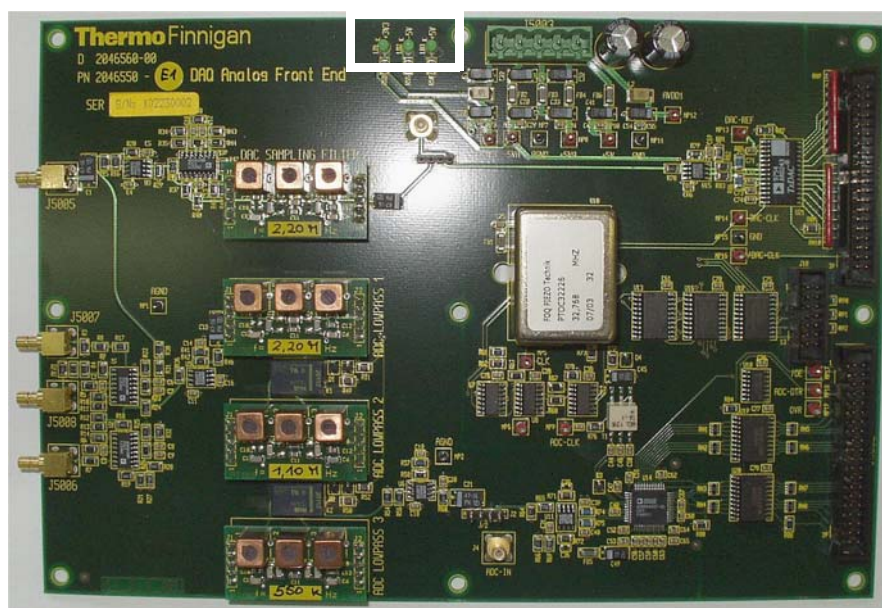


Figure 1-38. Data Acquisition Analog board

The diagnostic LEDs listed in Table 1-7 on page 1-51 show the status of the voltages applied to the board. The position of the LEDs on the board is indicated by a white rectangle in Figure 1-38.

Table 1-7. Diagnostic LEDs of the Data Acquisition Analog board

Name	Color	Description	Normal Operating Condition
+5 V	Green	+5 V voltage present	On
-5 V	Green	-5 V voltage present	On
+3.3 V	Green	+3.3 V voltage present	On

Power Supply 2 Board

The power supply 2 board provides the supply voltages for the data acquisition analog board. It is mounted to the back inside the housing of the internal computer. See Figure 1-36 on page 1-49.

The diagnostic LEDs listed in [Table 1-8](#) show the status of the voltages applied to the board.

Table 1-8. Diagnostic LEDs of the Power Supply 2 board

Name	Color	Description	Normal Operating Condition
+5.1 V	Green	+5.1 V voltage present	On
-5.1 V	Green	-5.1 V voltage present	On
+3.3 V	Green	+3.3 V voltage present	On

Instrument Control Board

[Figure 1-39](#) shows the instrument control board. The instrument control board is located in a housing next to the internal computer. It is connected to the Orbitrap Velos Pro MS main power.

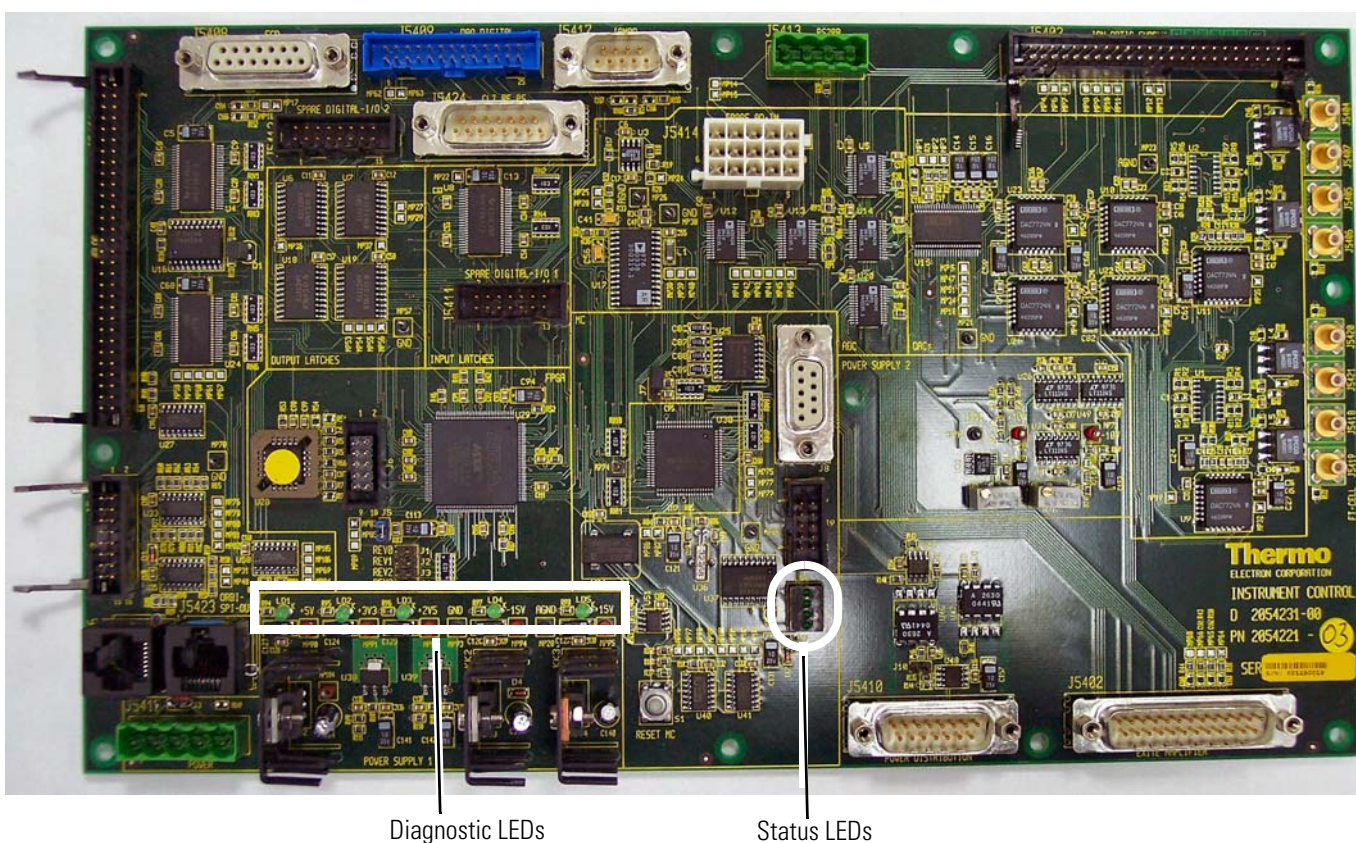


Figure 1-39. Instrument Control board

The instrument control board is used to interface the Velos Pro MS control electronics to the Orbitrap analyzer control electronics. Three signal lines are passed from the Velos Pro: a digital, parallel (DAC) bus, a serial SPI bus, and a Link Port Signal line. The instrument control board contains a micro controller, digital and analog converters, and serial port connectors.

On the instrument control board, analog signals from vacuum gauges are converted to digital signals and passed to the data system as well as to the power distribution board. (See [page 1-53](#).) Turbomolecular pumps (See “[Vacuum System](#)” on [page 1-30](#).) are attached to a serial port connector and this is connected via the signal lines to the linear ion trap.

The diagnostic LEDs listed in [Table 1-9](#) show the status of applied voltages to the board. The position of the diagnostic LEDs on the board is indicated by a white rectangle in [Figure 1-39](#) on [page 1-52](#).

Table 1-9. Diagnostic LEDs of the Instrument Control board

No.	Name	Color	Description	Normal Operating Condition
LD1	2.5 V	Green	2.55 V Input voltage present	On
LD2	3.3 V	Green	3.3 V Input voltage present	On
LD3	5 V	Green	5 V Input voltage present	On
LD4	-15 V	Green	-15 V Input voltage present	On
LD5	+15 V	Green	+15 V Input voltage present	On

Additionally, the board has four green LEDs that are directly connected to the micro controller. They indicate the state of the micro controller and possible error bits and can be used for software debugging. See [Table 1-10](#). The position of the status LEDs on the board is indicated by a white oval in [Figure 1-39](#) on [page 1-52](#).

Table 1-10. Software status LEDs of the Instrument Control board

No.	Description	Normal Operating Condition
6.1	Micro controller is working properly	Permanent flashing of LED
6.2	CAN bus connection to power distribution board enabled	On
6.3	Connection to internal computer and Velos Pro SPI bus enabled	On
6.4	Orbitrap analyzer SPI bus enabled	On Flashing on error

Power Distribution Board

[Figure 1-40](#) on [page 1-54](#) shows the power distribution board. It is located at the bottom of the right side of the instrument.

The power distribution board controls the vacuum system and the system power supplies, including the linear ion trap. Depending on the quality of the vacuum and the status of the turbomolecular pumps, it switches the vacuum gauges, the pumps, and the 230 V relays. It controls external relays with 24 V DC connections. In case of a vacuum failure, it initiates an automatic power down of the instrument.

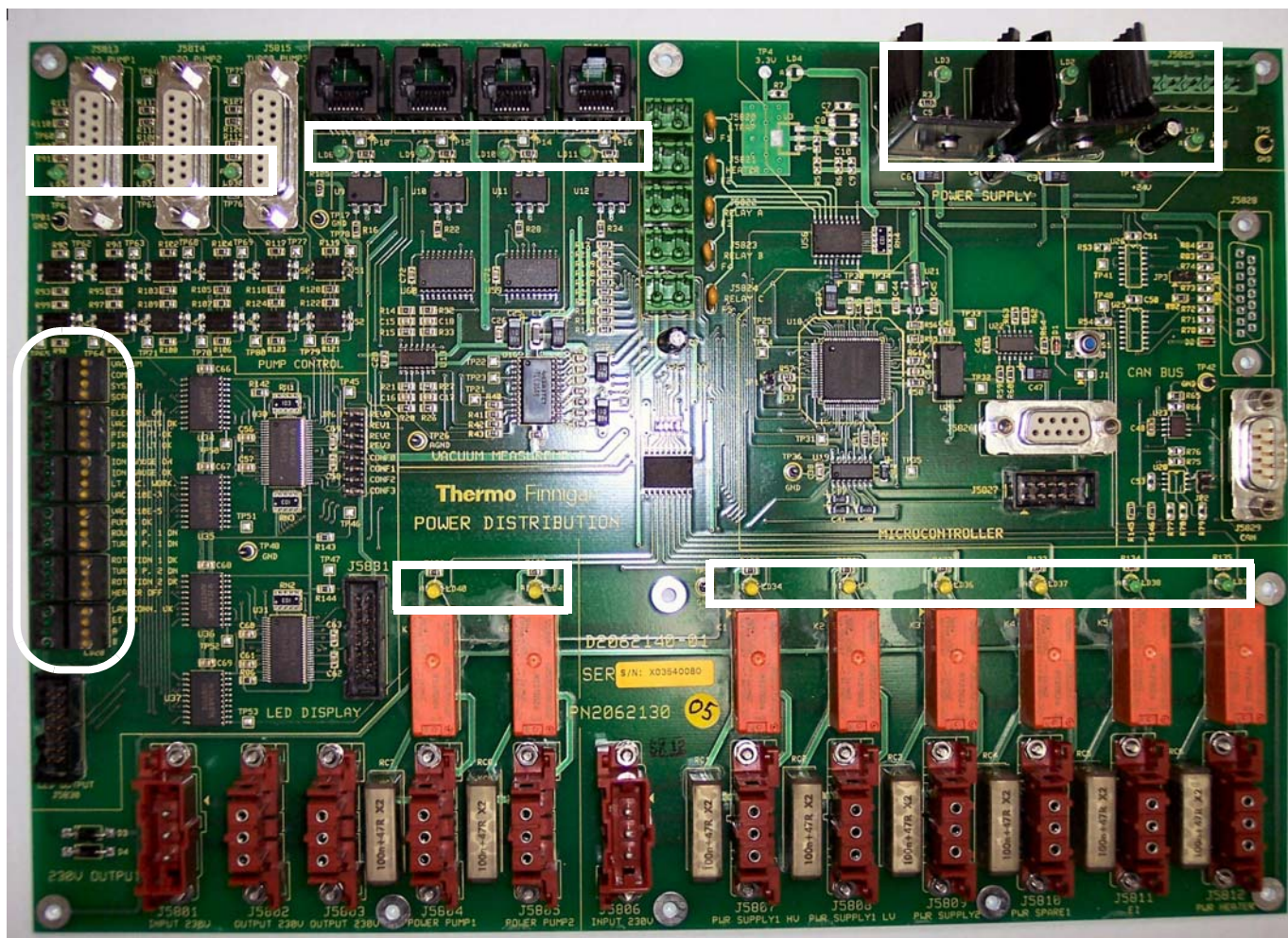


Figure 1-40. Power Distribution board

The power distribution board indicates all system states and error messages by status LEDs (See [Table 1-11](#) on [page 1-55.](#)) in the middle of the left side of the board. A green LED indicates that the status is OK. An orange LED indicates a status that differs from normal. The position of the LEDs on the board is indicated by a white oval in [Figure 1-40.](#)

The system status LEDs on the front side of the instrument (See [Figure 1-4](#) on [page 1-5.](#)) are controlled by the power distribution board. The information partially comes from external boards (for example, the Communication LED is controlled by the instrument control board). (See “[Instrument Control Board](#)” on [page 1-52.](#))

Diagnostic LEDs show the status of voltages applied from the board to other devices. The positions of the diagnostic LEDs on the board are indicated by white rectangles in [Figure 1-40.](#)

Table 1-11. Status LEDs of the Power Distribution board

LED green	LED orange	Information given by orange LED
Vacuum	High vacuum failure	High vacuum pressure > 10 ⁻⁸ mbar
Comm.	No communication with instrument control board	CAN bus problem or instrument control board not working
System	System is not ready	FT Electronics switch off or Vacuum Pumps switch off
Scan		Instrument is not scanning
Electr. On	Service mode	FT Electronics switch off
Vac. Units OK	Vacuum measurement failure	Vacuum gauge defective
Pirani Orbitrap analyzer OK	No function, at present	
Pirani LT OK	Pirani Velos Pro MS failure	Control signal < 0.5 V
Ion Gauge On	Penning Orbitrap Velos Pro MS Off	Forevacuum > 10 ⁻² mbar
Ion Gauge OK	Penning Orbitrap Velos Pro MS failure	Control signal < 0.5 V
LT Vacuum Work	Velos Pro MS vacuum failure	Vacuum forepump Velos Pro MS >10 ⁻¹ mbar
Vac. <10 ⁻³	Forevacuum failure	Forevacuum > 10 ⁻³ mbar
Vac. <10 ⁻⁵	High vacuum failure	High vacuum > 10 ⁻⁵ mbar
Pumps OK	Pumps Off	Pump down; leakage
Rough P. 1 On	Forepump #1 failure	Forepump defective
Turbo P. 1 On	TMP 1 failure	TMP defective/error ^a
Rotation 1 OK	TMP 1 failure	80% rotation speed of TMP not reached
Turbo P. 2 On	TMP 2 failure	TMP defective/error ^a
Rotation 2 OK	TMP 2 failure	80% rotation speed of TMP not reached
Heater Off	Heater enabled	Heater enabled
LAN Conn. OK	LAN connection failure	LAN interrupted (Option)
EI On	No function, at present	
A	System reset	System reset has occurred
B		Micro controller idle

^aAn error of TMP 3 is indicated by an LED directly located on the pump controller. An error of TMP 4 is indicated in the software.

Depending on user actions, the power distribution is switched to various working modes by the hardware. See [Table 1-12](#) on [page 1-56](#).

Table 1-12. Working modes of the Power Distribution board

Action	Consequences
a. Main switch off	Complete system including linear ion trap and multiple socket outlets (ETD Module, for example) are without power
b. Vacuum Pumps switch off	Everything is switched off
c. FT Electronics switch off	All components are switched off with exception of the following ones: <ul style="list-style-type: none"> • Heater control • Multiple socket outlets • Power distribution board • Pumps • Vacuum control • Velos Pro MS (has a separate Service switch)

Table 1-13 shows the possible operating states of the power distribution.

Table 1-13. Operating states of the Power Distribution board

Action	Consequences
1. Main switch on, Vacuum Pumps switch off	Everything is switched off
2. Vacuum Pumps switch on and FT Electronics switch on	System starts up: pumps and electronics switched on
3. Check linear ion trap and Orbitrap Velos Pro MS forevacuum pumps: 10^{-0} mbar after 30 s.	If not ok: switch off system and light error LED [®] ; power distribution remains switched on
4. After the system has started, the Pirani gauge returns a vacuum $< 10^{-2}$ mbar and both TMPs reach 80% rotation speed	Switch on Penning gauge
5. Vacuum and 80% rotation speed of TMPs not reached after preset time (< 8 min, otherwise the pumps automatically switch off).	Switch off system (including linear ion trap) and light error LED [®] ; power distribution remains switched on
6. One or more vacuum gauges defective (control signal < 0.5 V).	Light error LED only, otherwise ignore
7. After the operating status is reached, the pressure at one gauge exceeds the security threshold for more than the preset time period: <ul style="list-style-type: none"> • Pirani gauge Orbitrap Velos Pro MS $> 10^{-1}$ mbar • Penning gauge Orbitrap Velos Pro MS $> 10^{-3}$ mbar • Pirani gauge Velos Pro MS forepump $> 10^{-1}$ mbar 	System is shut down with exception of power distribution (light error LED). Rebooting of the system by switching off/on of the main switch.

Table 1-13. Operating states of the Power Distribution board, continued

Action		Consequences
8.	Rotation speed of a TMP falls below 80%	Shut down system (see 7.); light LED* of corresponding pump.
9.	Service switch linear ion trap off	Linear ion trap electronics switched off, pumps keep on running; Orbitrap Velos Pro MS without data link, keeps on running
10.	FT Electronics switch Orbitrap Velos Pro MS off	Orbitrap Velos Pro MS electronics switched off, pumps keep on running; Orbitrap Velos Pro MS without data link, keeps on running
11.	Failure of linear ion trap or Orbitrap Velos Pro MS (for example, fuse is opened).	If the vacuum in one part deteriorates, the complete system is shut down.
12.	Mains failure	System powers up after the electricity is available again. All devices reach the defined state. Linear ion trap and internal computer must reboot.

*After the shutdown, the LED flashes that represents the reason for the shutdown.

Power Supply 1 Board

Figure 1-41 shows the power supply 1 board. This board is located next to the power distribution board. It provides the power for the ion optic supply board (See “Ion Optic Supply Board” on page 1-60.) and the instrument control board. (See “Instrument Control Board” on page 1-52.)

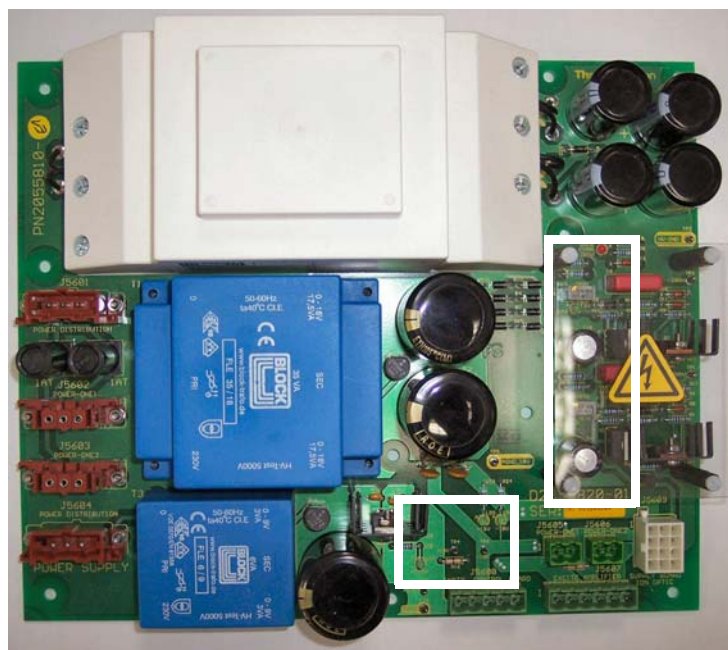


Figure 1-41. Power Supply 1 board



Warning Electrical Shock Hazard. Parts of the power supply 1 board are at high voltage. Shut down the instrument and disconnect it from line power before performing service. ▲

The diagnostic LEDs listed in Table 1-14 show the status of the voltages applied to the board. The position of the LEDs on the board is indicated by the white rectangles in Figure 1-41.

Table 1-14. Diagnostic LEDs of the Power Supply 1 board

Name	Color	Description	Normal Operating Condition
+285 V	Green	+285 V Output voltage present	On
-285 V	Green	-285 V Output voltage present	On
Over Current +285 V	Red	LED lit dark red: $I_{out} > 80$ mA LED lit bright red: output is short-circuited	Off
Over Current -285 V	Red	LED lit dark red: $I_{out} > 80$ mA LED lit bright red: output is short-circuited	Off

Table 1-14. Diagnostic LEDs of the Power Supply 1 board, continued

Name	Color	Description	Normal Operating Condition
+18 V	Green	+18 V Output voltage present	On
-18 V	Green	-18 V Output voltage present	On
+8.5 V	Green	+8.5 V Output voltage present	On

Electronic Boards on the Left Side of the Instrument

Figure 1-42 shows the left side of the instrument with the panel opened. This side of the instrument contains mostly boards that are part of the Orbitrap analyzer control.

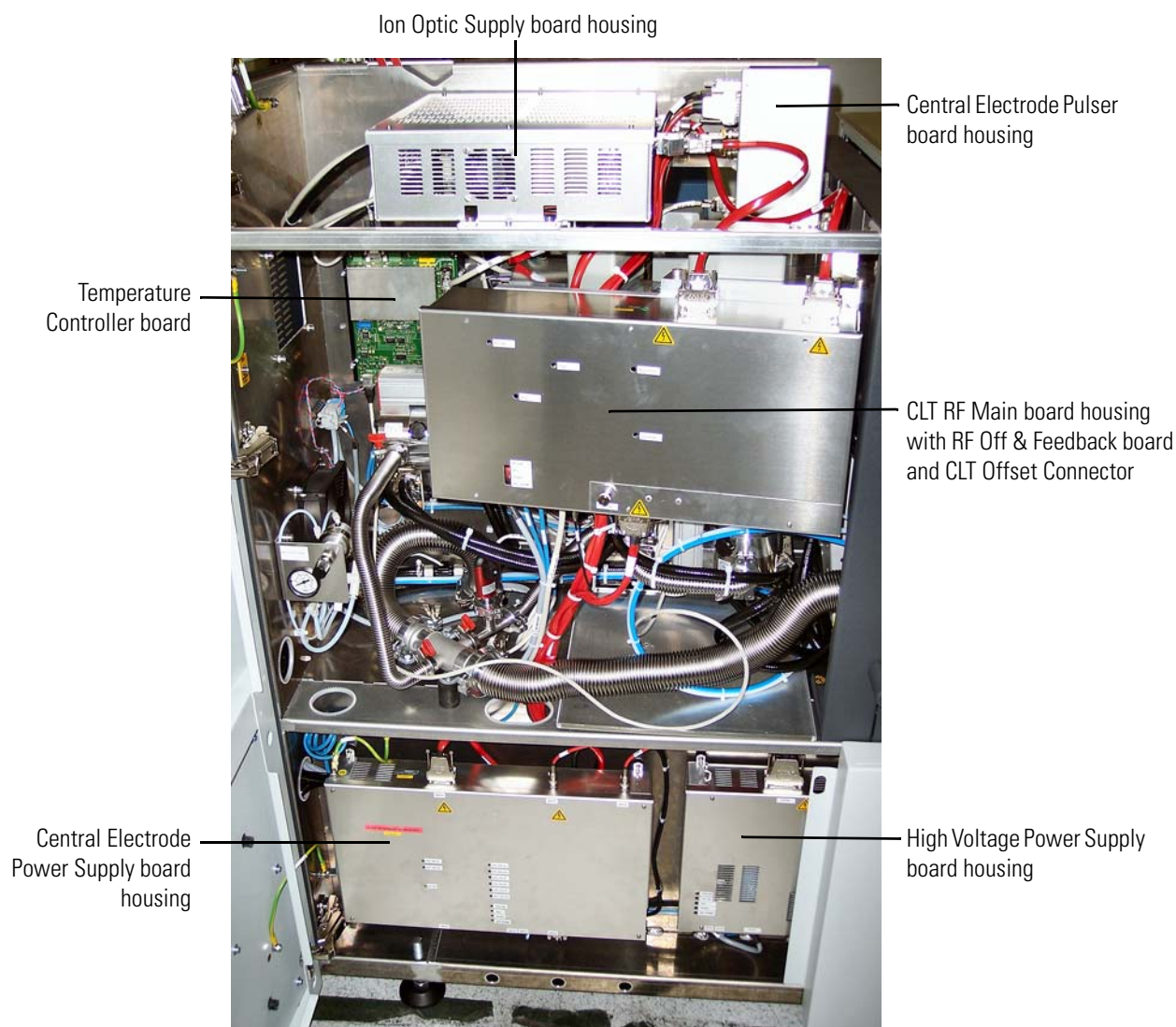


Figure 1-42. Electronic boards on the left side of the instrument

The main components on this side are described starting from the top.

Ion Optic Supply Board

Figure 1-43 on page 1-60 shows the ion optic supply board. The board is located in a housing on top of the left instrument side of the instrument. This board supplies the voltages and the radio frequency for the ion guides and interoctapole lenses of the Orbitrap Velos Pro mass spectrometer. It has an RF detector for the RF output control. The board also provides the entrance voltage, the exit voltage, and the reflector DC voltages as well as the RF voltages to the octapole of the Orbitrap analyzer. See “Orbitrap Analyzer” on page 1-14 for further information.



Figure 1-43. Ion Optic Supply board

The diagnostic LEDs listed in Table 1-15 show the status of applied voltages to the board. The position of the LEDs on the board is indicated by white rectangles in Figure 1-43.



Warning Electrical Shock Hazard. Parts of the board are at high voltage. Shut down the instrument and disconnect it from line power before performing service. ▲

Table 1-15. Diagnostic LEDs of the Ion Optic Supply board

No.	Name	Color	Description	Normal Operating Condition
LD1	+275 V	Green	+275 V Input voltage present	On
LD2	-275 V	Green	-275 V Input voltage present	On

Table 1-15. Diagnostic LEDs of the Ion Optic Supply board, continued

No.	Name	Color	Description	Normal Operating Condition
LD3	+29 V	Green	+29 V Input voltage present	On
LD5	+15 V	Green	+15 V Input voltage present	On
LD6	-15 V	Green	-15 V Input voltage present	On
LD7	RF1_ON	Blue	RF1 generator switched on	depending on application; LED flashes during scanning
LD8	RF2_ON	Blue	RF2 generator switched on	depending on application; LED flashes during scanning

Central Electrode Pulser Board

The central electrode pulser board is located in a housing that is mounted to the flange of the UHV chamber. See [Figure 1-44](#).

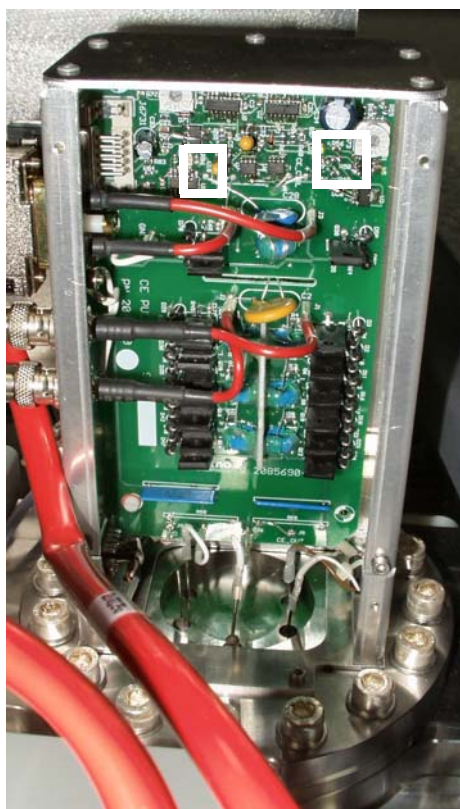


Figure 1-44. Central Electrode Pulser board

The board switches the injection and measurement voltages for the central electrode and the detection electrodes of the Orbitrap analyzer. Resistor-capacitor circuits on the board convert the switching pulse into a smooth transition between the voltages.

The diagnostic LEDs listed in [Table 1-16](#) on [page 1-62](#) show the status of the voltages applied to the board as well as some operating states. The position of the LEDs on the board is indicated by the white rectangles in [Figure 1-44](#).

Table 1-16. Diagnostic LEDs of the Central Electrode Pulser board

No.	Name	Color	Description	Normal Operating Condition
LD1	TRIG	Green	Trigger signal indicator	Flashing when scanning
LD2	PS	Green	24V Power Supply is OK	On

Temperature Controller Board

The temperature controller board is located on the top left side of the instrument, next to the CLT RF main board. See [Figure 1-42](#) on [page 1-59](#). The temperature controller board keeps the temperature of the analyzer chamber to a preset value. A Peltier element that can be used for heating as well as for cooling is used as an actuator. Activation is done via the serial SPI (Serial Peripheral Interface) bus.

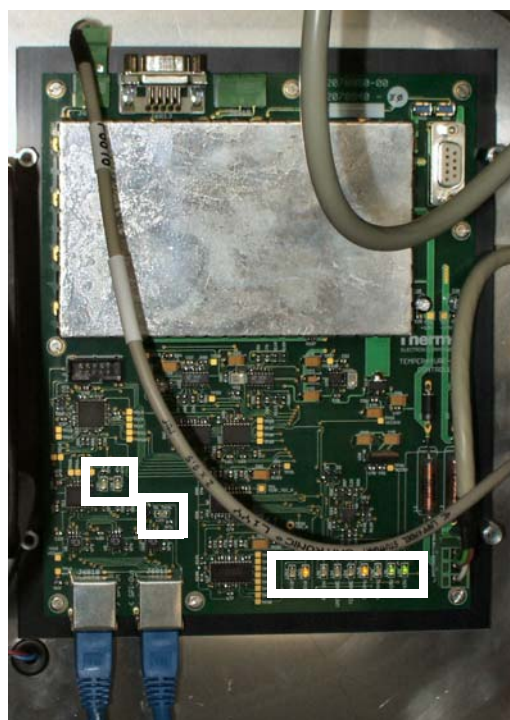


Figure 1-45. Temperature Controller board

The diagnostic LEDs listed in [Table 1-17](#) on [page 1-63](#) show the status of the voltages applied to the board as well as some operating states. The positions of the LEDs on the board are indicated by the white rectangles in [Figure 1-45](#).

Table 1-17. Diagnostic LEDs of the Temperature Controller board

No.	Name	Color	Description	Normal Operating Condition
LD1	+15 V	Green	+15 V Input voltage present	On
LD2	-15 V	Green	-15 V Input voltage present	On
LD3	TEC >60C	Yellow	Temperature of cold side Peltier element above 60 °C	Off
LD4	Unit >60C	Yellow	Temperature of UNIT heat sink above 60 °C	Off
LD5	Reg Off	Yellow	Control switched off	Off
LD6	No Term	Yellow	SPI bus termination board missing	Off
LD7	SDT enable	Green	Interface has been addressed and sends/receives data	Flashing on SPI bus data transfer
LD8	SEL	Green	Board has been addressed	Flashing on SPI bus data transfer
LD9	Heating	Yellow	Peltier element is heating	Depending on system state
LD10	Cooling	Yellow	Peltier element is cooling	Depending on system state
LD11	UR>0	Yellow	Summation voltage controller >0 V	Off when adjusted
LD12	UR<0	Yellow	Summation voltage controller <0 V	Off when adjusted

CLT RF Unit

The CLT RF unit comprises the CLT RF main board and the RF off & feedback board. The unit operates the curved linear trap (CLT) with four phases RF voltage and three pulsed DC voltages (PUSH, PULL, and OFFSET).

The CLT RF main board is located in a housing in the center of the left side of the instrument. See [Figure 1-42](#) on [page 1-59](#). This board provides an RF voltage (“Main RF”) for the curved linear trap. It allows switching off the RF and simultaneous pulsing of each CLT electrode. See “[Orbitrap Analyzer](#)” on [page 1-14](#) for further information. The board communicates with the instrument control board via an SPI bus.

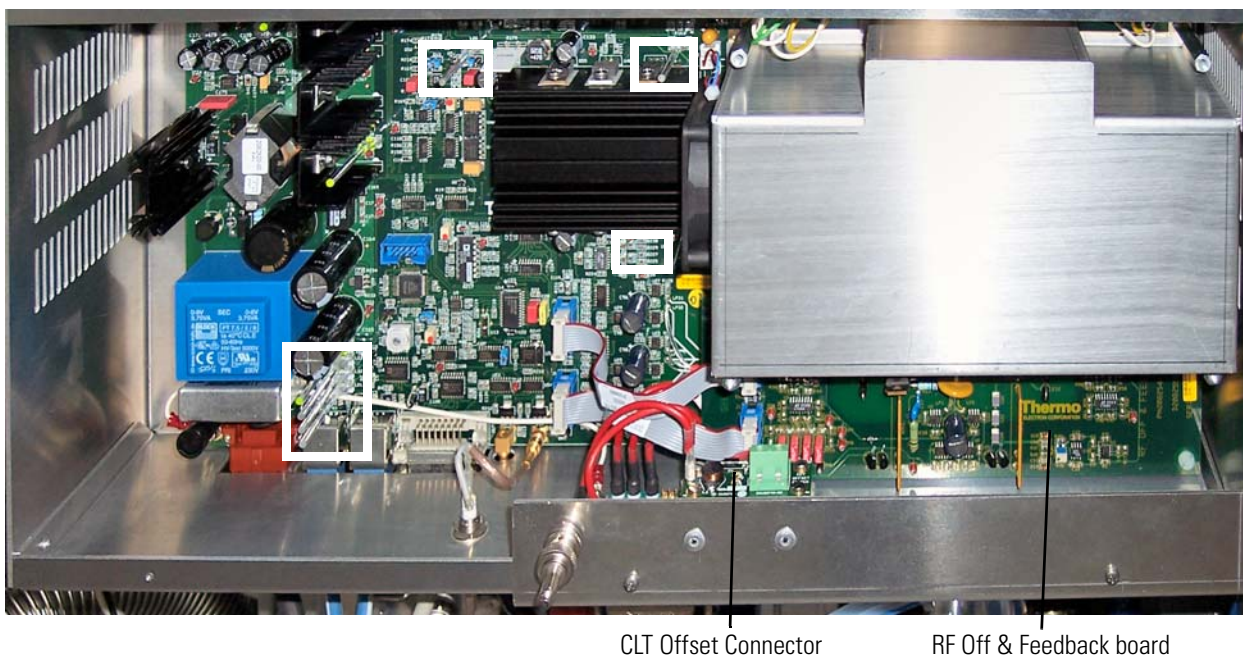


Figure 1-46. CLT RF unit (cover removed)

The diagnostic LEDs listed in [Table 1-18](#) show the status of the voltages applied to the board as well as some operating states. The position of the LEDs on the board is indicated by the white rectangles in [Figure 1-46](#).

Table 1-18. Diagnostic LEDs of the CLT RF Main board

No.	Name	Color	Description	Normal Operating Condition
LD1	NO TERM	Yellow	SPI bus termination board missing	Off
LD2	SEND	Yellow	Interface has been addressed and sends/receives data	Flashing on SPI bus data transfer
LD3	SEL	Green	Board has been addressed	Flashing on SPI bus data transfer
LD4	RF ON	Green	RF voltage on	On
LD5	NO LOCK	Yellow	PLL has been not locked	50% intensity
LD6	OVL	Yellow	RF Amplifier overload	Off
LD7	OVHEAT	Red	Heatsink temperature > 73 °C	Off

The RF off & feedback board is an add-on board to the CLT RF main board. It is located in the same housing. See [Figure 1-46](#) on [page 1-64](#).

The CLT Offset connector, which removes interfering signals from the circuit, is also mounted in the housing.

Central Electrode Power Supply Board

The central electrode power supply board is mounted in a housing on the bottom left side of the instrument. See [Figure 1-47](#).

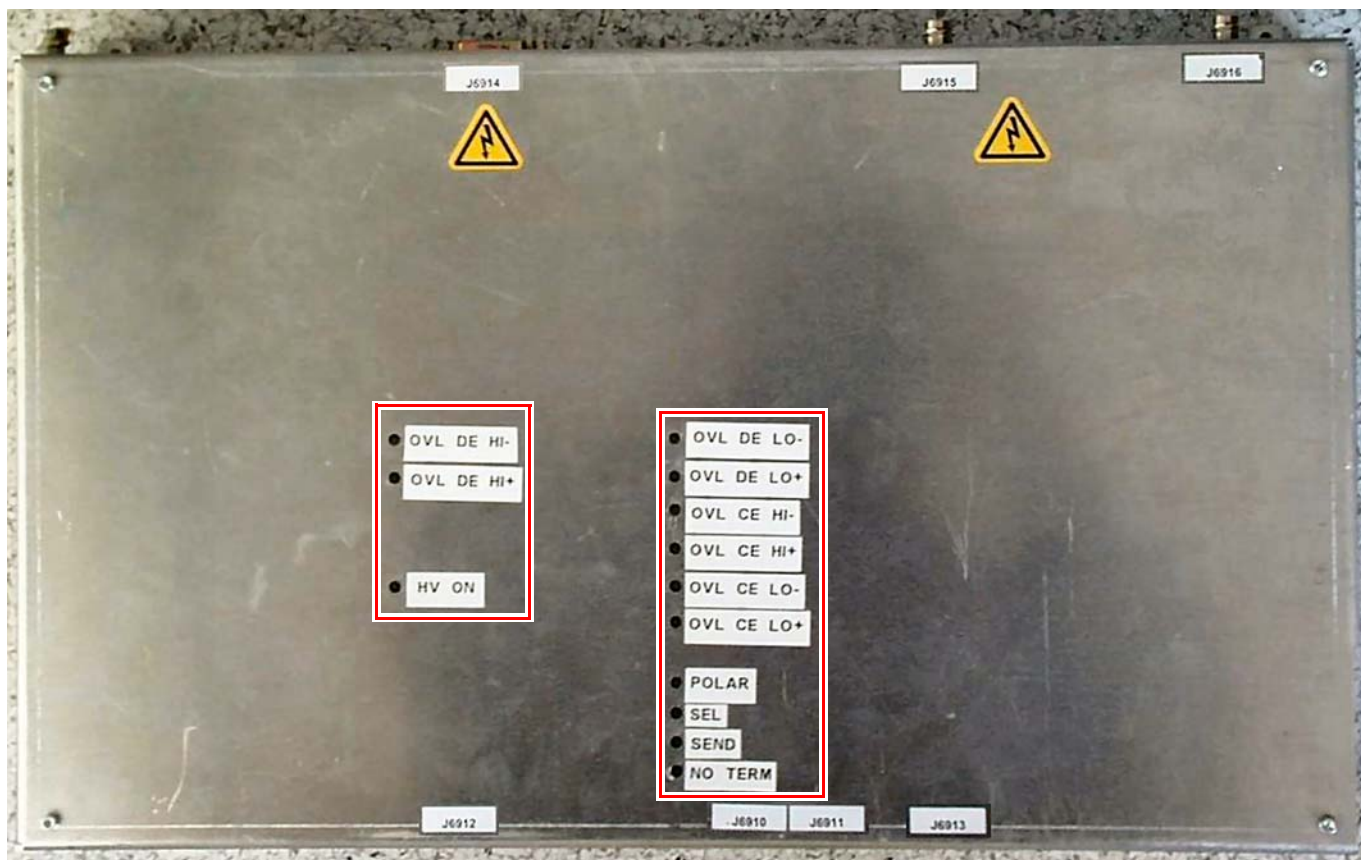


Figure 1-47. Central Electrode Power Supply board

The board supplies four DC voltages to the Orbitrap analyzer:

- Two central electrode (CE) voltages: CE HIGH and CE LOW.
- Two deflector electrode (DE) voltages: DE HIGH and DE LOW.

For positive ions, the CE voltages are negative and the DE voltages are positive. The maximum CE voltage is 3 kV and the maximum DE voltage is 1 kV. The board communicates via the SPI bus.

In addition to a ventilator on the bottom right side, a water-cooled Peltier element on the rear side of the board serves as means of heat dissipation.

The diagnostic LEDs listed in [Table 1-19](#) show the status of the voltages applied to the board as well as some operating states. The position of the LEDs on the board is indicated by the red rectangles in [Figure 1-47](#) on [page 1-65](#).

Table 1-19. Diagnostic LEDs of the Central Electrode Power Supply board

No.	Name	Color	Description	Normal Operating Condition
LD1	OVL DE HI-	Yellow	Negative side of Deflector High Supply has been overloaded	Off when HV is switched on
LD2	OVL DE HI+	Yellow	Positive side of Deflector High Supply has been overloaded	Off when HV is switched on
LD3	No Term	Red	SPI bus termination board missing	Off
LD4	Send	Yellow	Interface has been addressed and sends/receives data	Flashing on SPI bus data transfer
LD5	Sel	Green	Board has been addressed	Flashing on SPI bus data transfer
LD6	Polarity	Blue	Positive/negative ion mode	Off (positive mode)
LD7	OVL CE LO+	Yellow	Positive side of Central Electrode Low Supply has been overloaded	Off when HV is switched on
LD8	OVL CE LO-	Yellow	Negative side of Central Electrode Low Supply has been overloaded	Off when HV is switched on
LD9	OVL CE HI+	Yellow	Positive side of Central Electrode High Supply has been overloaded	Off when HV is switched on
LD10	OVL CE HI-	Yellow	Negative side of Central Electrode High Supply has been overloaded	Off when HV is switched on
LD11	OVL DE LO+	Yellow	Positive side of Deflector Low Supply has been overloaded	Off when HV is switched on
LD12	OVL DE LO-	Yellow	Negative side of Deflector Low Supply has been overloaded	Off when HV is switched on
LD13	HV ON	Green	High voltage switched on	On when HV is switched on

High Voltage Power Supply Board

The high voltage power supply board is mounted in a housing on the bottom left side of the instrument. See [Figure 1-42](#) on [page 1-59](#). This board provides five DC voltages for the ion optics of the Orbitrap Velos Pro mass spectrometer. Two voltages supply the lenses of the instrument. Three voltages are applied to the RF CLT main board to be

used as focusing potentials for the curved linear trap. See “Orbitrap Analyzer” on page 1-14 for further information. The board communicates via the SPI bus.



Warning Electrical Shock Hazard. The high voltage power supply board creates voltages up to 3.5 kV! Shut down the instrument and disconnect it from line power before performing service. ▲

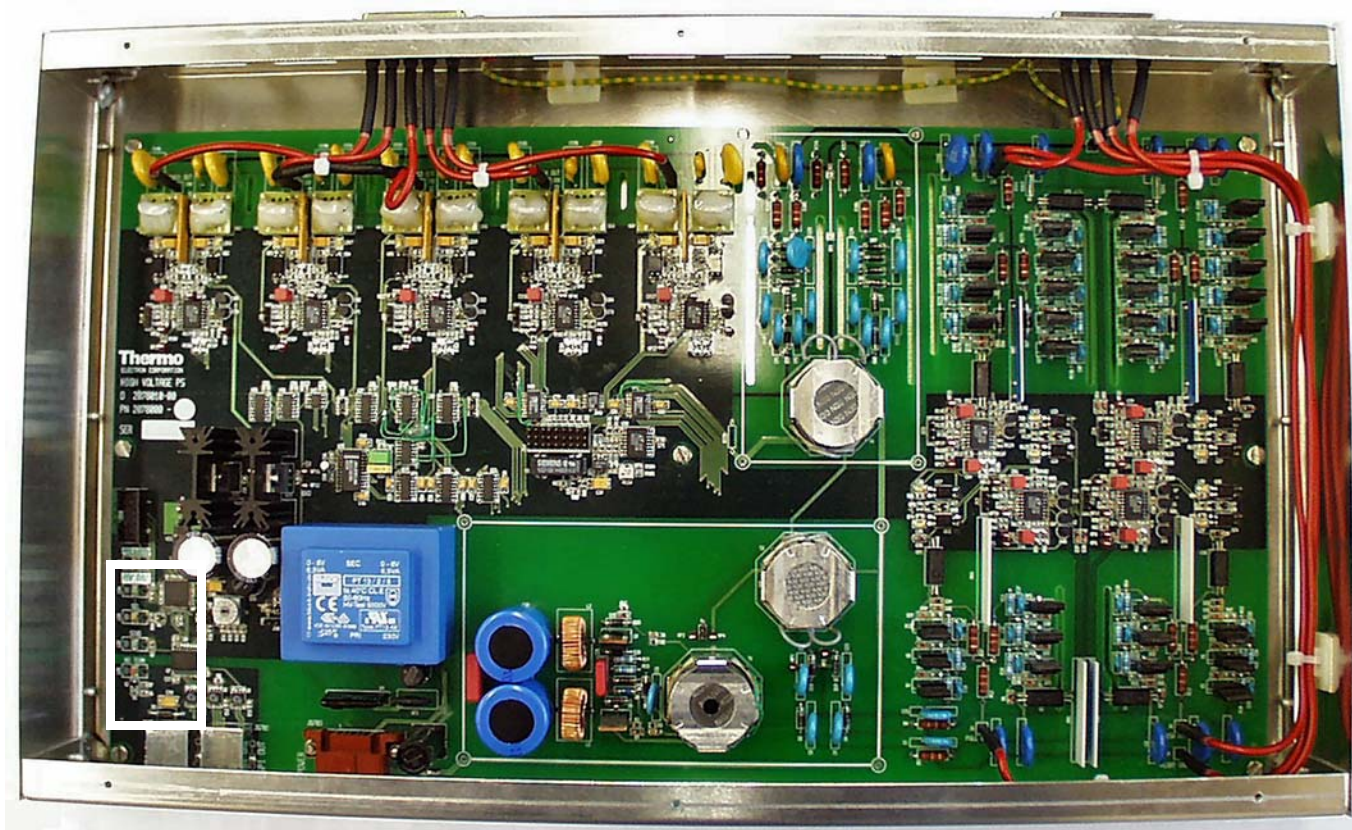


Figure 1-48. High Voltage Power Supply board (cover removed)

The diagnostic LEDs listed in Table 1-20 on page 1-68 show the operating states of the board. The position of the LEDs on the board is indicated by the white rectangles in Figure 1-48.

Table 1-20. Diagnostic LEDs of the High Voltage Power Supply board

No.	Name	Color	Description	Normal Operating Condition
LD1	NO TERM	Red	SPI bus termination board missing	Off
LD2	SEND	Yellow	Interface has been addressed and sends/receives data	Flashing on SPI bus data transfer
LD3	SEL	Green	Board has been addressed	Flashing on SPI bus data transfer
LD4	HV ON	Green	High voltage is switched on	On
LD5	POLARITY	Green	Positive/negative ion mode	Off (positive mode)

SPI Bus Termination Board

Various boards communicate via the SPI bus, a serial RS485-based bus system. The SPI Bus Termination board reduces unwanted signal reflections. The boards indicate a missing termination (after maintenance, for example) by LEDs.

The SPI Bus Termination board is located below the High Voltage Power Supply board, at the bottom left side of the instrument. See [Figure 1-49](#).



SPI bus termination board

Figure 1-49. High Voltage Power Supply board with SPI Bus Termination board

Chapter 2 Basic System Operations

Many maintenance procedures for the Orbitrap Velos Pro system require that the mass spectrometer be shut down. In addition, the Orbitrap Velos Pro system can be placed in Standby condition if the system is not to be used for 12 hours or more.

The following topics are described in this chapter:

- “Shutting Down the System in an Emergency” on page 2-2
- “Placing the Instrument in Standby Condition” on page 2-4
- “Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely” on page 2-7
- “Starting Up the System after a Shutdown” on page 2-9
- “Resetting the System” on page 2-12
- “Resetting Tune and Calibration Parameters to their Default Values” on page 2-13
- “Turning Off the Reagent Ion Source: What to Expect” on page 2-14

Shutting Down the System in an Emergency

If you need to turn off the mass spectrometer in an emergency, place the main power switch (located on the power panel at the right side of the Orbitrap Velos Pro mass spectrometer) in the Off (0) position. This turns off all power to the instrument, including the linear ion trap, multiple socket outlets, and the vacuum pumps. The main power switch must be turned 90° anti-clockwise to switch off the instrument. See [Figure 2-1](#).

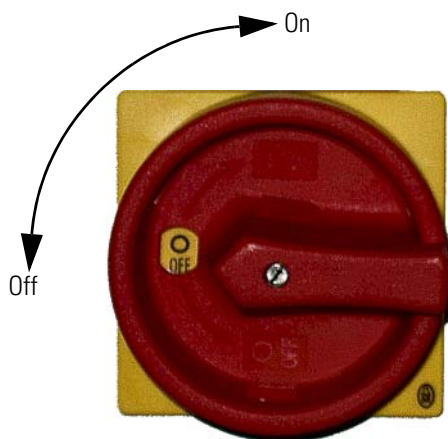


Figure 2-1. Main power switch in Off position

The instrument is automatically vented by the vent valve of the linear ion trap. The vent valve vents the system 30 seconds after power is switched off.

Although removing power abruptly will not harm any component within the system, this is not the recommended shutdown procedure to follow. See [“Shutting Down the Instrument”](#) on [page 2-7](#) for the recommended procedure.

Note To separately turn off the recirculating chiller or computer in an emergency, use the On/Off switches on the chiller and computer, respectively. ▲

Behavior of the System in Case of a Main Failure

A main power failure has the same consequence as switching off via the main power switch. If the power is available again, the system starts up automatically: the pumps are switched on and the instrument is pumped down. If the system has been vented during the mains failure, it is necessary to bake out the system to obtain the operating vacuum. See [“Baking Out the System”](#) on [page 3-4](#).

It is not possible to check whether the system was vented. The log file of the data system indicates a reboot of the system. In case of frequent but short power failures we recommend installing an uninterruptible power supply (UPS). If main power failures occur frequently while the system is not attended (for example, in the night), we recommend installing a power fail detector.

Note The intentional venting of the system is performed with the vent valve of the linear ion trap. ▲

Placing the Instrument in Standby Condition

The Orbitrap Velos Pro system should not be shut down completely if you are not going to use it for a short period of time, such as overnight or over the weekend. When you are not going to operate the system for 12 hours or more, you can leave the system in Standby condition.

In case of an Orbitrap Velos Pro ETD mass spectrometer, first place the ETD Module in Standby condition. Then place the mass spectrometer in Standby condition according to the procedures described in the following topics.

Placing the ETD Module in Standby Condition

❖ To place the ETD Module in Standby condition

1. If the Tune Plus window is not already open, choose **Start > Programs > Thermo Instruments > LTQ > LTQ Tune** from the taskbar. The Tune Plus window will open.

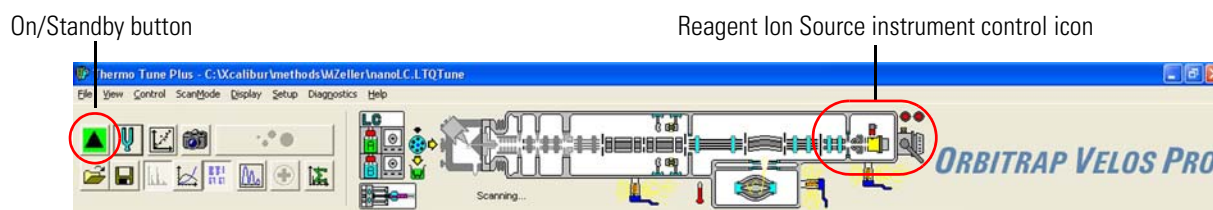


Figure 2-2. Tune Plus window (Orbitrap Velos Pro ETD), toolbar



You can determine the state of the MS detector by observing the state of the On/Off/Standby button on the Control/Scan Mode toolbar. See [Figure 2-2](#). The three different states of the On/Standby button are shown at the left.

2. Click the Reagent Ion Source portion of the instrument control graphic at the top of the Tune Plus window. (See [Figure 2-2](#).) The Reagent Ion Source dialog box appears. (See [Figure 2-3](#).)
3. In the Reagent Ion Source dialog box, clear the Reagent Ion Source On box to place the Reagent Ion Source in Standby condition. See [Figure 2-3](#) on [page 2-5](#). This places the Reagent Ion Source in Standby condition as indicated by the Actual condition shown to the right of the Reagent Ion Source On box.

When the reagent ion source is placed in Standby condition, the filament and vial heaters turn off. Simultaneously, a valve opens to allow the nitrogen gas to cool the reagent vials. This cooling nitrogen runs until the reagent vials reach 70 °C. The audible rush (hissing noise) of nitrogen from the reagent ion source area in the back of the ETD Module is normal operation.

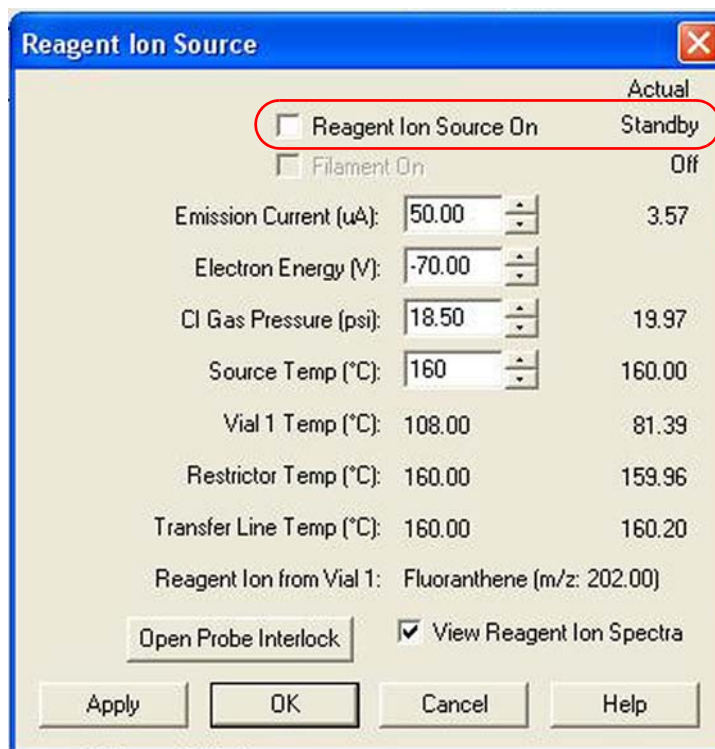


Figure 2-3. Reagent Ion Source dialog box with Reagent Ion Source On box and Actual condition circled

If the reagent ion source is on when you place the Orbitrap Velos Pro ETD mass spectrometer in Standby mode, the filament turns off immediately. In contrast, the vial heaters stay on for 60 minutes before they turn off and the cooling gas begins. Because the filament is turned off, you can perform minor maintenance procedures on the ETD Module without cooling the reagent inlet.



Warning Burn Hazard. Install or exchange the reagent vials by following the procedure in “Replacing the Reagent Vials” on page 3-48. The reagent vials will be too hot to touch after the cooling nitrogen turns off at 70°C. Verify that the reagent vials are cool to the touch before handling them. ▲

More information about turning on and off the reagent heaters is given in “Reagent Heaters” on page 1-27.



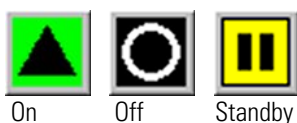
Warning Burn Hazard. The restrictor, the transfer line, and the ion source heater operate at 160 °C. Do not attempt to touch them unless the Orbitrap Velos Pro mass spectrometer is shut down (See “Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely” on page 2-7.) and these heaters have had sufficient time to cool down to room temperature. ▲

Placing the MS in Standby Condition

❖ **To place the Orbitrap Velos Pro system in Standby condition**

1. Wait until data acquisition, if any, is complete.
2. Turn off the flow of solvent from the LC (or other sample introduction device).

Note For instructions on how to operate the LC from the front panel, refer to the manual that came with the LC. ▲



3. From the Tune Plus window, choose **Control > Standby** (or click the **On/Standby** button to toggle it to Standby) to put the instrument in Standby condition. The consequences of this user action are described in the *LTQ Series Hardware Manual*. The System LED on the front panel of the LTQ Velos Pro mass spectrometer turns yellow when the system is in Standby condition.
4. Leave the LC power on.
5. Leave the autosampler power on.
6. Leave the data system power on.
7. Leave the Orbitrap Velos Pro MS main power switch in the On position.

Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely

The Orbitrap Velos Pro mass spectrometer does not need to be shut down completely if you are not going to use it for a short period of time, such as overnight or over weekends. Shut down ETD Module and MS system completely only if you do not want to use them for an extended period or if you want to perform a maintenance or service procedure.

❖ To shut down the instrument completely

1. Place the ETD Module in Standby condition as described in [“Placing the ETD Module in Standby Condition”](#) on [page 2-4](#).
2. Shut down the instrument as described in [“Shutting Down the Instrument”](#) below. This also shuts down the ETD Module because its power controls are linked to the Orbitrap Velos Pro MS power controls through the ETD Module Interface board. See [“ETD Module Interface Board”](#) on [page 1-25](#).

Shutting Down the Instrument

❖ To shut down the Orbitrap Velos Pro system

1. Wait until data acquisition, if any, is complete.
2. Turn off the flow of solvent from the LC (or other sample introduction device).

Note For instructions on how to operate the LC from the front panel, refer to the manual that came with the LC. ▲

3. From the Tune Plus window, choose **Control > Off** to put the instrument in Off condition. When you choose **Control > Off**, all high voltages are shut off, as are the flows of the sheath gas and the auxiliary gas.
4. Put the FT Electronics switch to the Off position. See [Figure 1-7](#) on [page 1-8](#).
5. Put the Vacuum Pumps switch to the Off position. See [Figure 1-7](#). When you place the switch in the Off position, the following occurs:
 - a. All power to the instrument, including the turbomolecular pumps and the rotary-vane pumps, is turned off.
 - b. After 30 s, power to the vent valve solenoid of the ion trap is shut off. The vent valve opens and the vacuum manifold is vented with nitrogen to atmospheric pressure through a filter. You can hear a hissing sound as the gas passes through the filter.

Basic System Operations

Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely

6. Leave the main power switch of the Orbitrap Velos Pro mass spectrometer in the On position.
7. During service or maintenance operations that require opening the vacuum system of the LTQ Velos Pro MS or the Orbitrap Velos Pro MS, always put the main switch (main circuit breaker) to the Off position. You can secure the main switch with a padlock or tie-wrap to prevent unintended re-powering.



Warning Burn Hazard. Allow heated components to cool down before you service them (the ion transfer tube is operated at about 300 °C, for example). ▲

Note If you are planning to perform routine or preventive system maintenance on the Orbitrap Velos Pro mass spectrometer only, you do not need to turn off the recirculating chiller, LC, autosampler, or data system. In this case, the shutdown procedure is completed. However, if you do not plan to operate your system for an extended period of time, you might want to turn off the recirculating chiller, LC, autosampler, and data system. ▲

Starting Up the System after a Shutdown

To start up the Orbitrap Velos Pro mass spectrometer after it has been shut down, you need to do the following:

1. Start up the instrument.
2. Set up conditions for operation.

Starting Up the Instrument

Note The recirculating chiller and data system must be running before you start up the instrument. The instrument will not operate until it has established a communication link to the data system. ▲

❖ To start up the Orbitrap Velos Pro mass spectrometer

1. Start up the (optional) LC and autosampler as is described in the manual that came with the LC and autosampler.
2. Start up the data system and the chiller.
3. Turn on the flows of helium and nitrogen at the tanks, if they are off.
4. Make sure that the main power switch of the LTQ Velos Pro MS is in the On position and the electronics service switch of the LTQ Velos Pro MS is in the Operating position.
5. Place the main power switch at the right side of the Orbitrap Velos Pro mass spectrometer in the On position.
6. Put the Vacuum Pumps switch to the On position. See [Figure 1-7](#) on [page 1-8](#). The rotary-vane pumps and the turbomolecular pumps are started.

Note Pumping the system after a complete shut down takes hours and requires overnight baking of the system. ▲

7. Put the FT Electronics switch to the On position. See [Figure 1-7](#). When you place the FT Electronics switch to the On position, the following occurs:
 - a. Power is provided to all electronic boards. (The electron multiplier, conversion dynode, 8 kV power to the API source, main RF voltage, and quadrupole RF voltage remain off.)
 - b. The internal computer reboots. After several seconds, the Communication LED on the front panel turns yellow to

indicate that the data system has started to establish a communication link.

- c. After several more seconds, the Communication LED turns green to indicate that the data system has established a communication link. Software for the operation of the instrument is then transferred from the data system to the instrument.
- d. After three minutes, the System LED of the ion trap turns yellow to indicate that the software transfer from the data system is complete and that the instrument is in Standby condition.

Note The Vacuum LED on the front panel of the LTQ Velos Pro turns green only if the pressure in the vacuum manifold is below the maximum allowable pressure (5×10^{-4} Torr in the analyzer region, and 2 Torr in the S-lens region), and the safety interlock switch on the API source is pressed down (that is, the API flange is secured to the spray shield). ▲

8. Press the Reset button on the LTQ Velos Pro to establish the communication link between LTQ Velos Pro and internal computer.

If you have an LC or autosampler, start it as is described in the manual that came with the LC or autosampler. Then, proceed to “[Setting Up Conditions for Operation](#)“. If you do not have either, go to the topic directly.

Setting Up Conditions for Operation

❖ To set up your Orbitrap Velos Pro mass spectrometer for operation

1. Before you begin data acquisition with your Orbitrap Velos Pro system, you need to allow the system to pump down for at least eight hours. Operation of the system with excessive air and water in the vacuum manifold can cause reduced sensitivity, tuning problems, and a reduced lifetime of the electron multiplier.

Note The vacuum in the analyzer system can be improved by an overnight baking of the system. See “[Baking Out the System](#)” on [page 3-4](#). ▲

2. Ensure that the gas pressures are within the operational limits:
 - Helium: 275 ± 70 kPa (2.75 ± 0.7 bar, 40 ± 10 psi),
 - Nitrogen: 690 ± 140 kPa (6.9 ± 1.4 bar, 100 ± 20 psi).

In case of an Orbitrap Velos Pro ETD mass spectrometer, also check the pressure of the reagent carrier gas: 690 ± 140 kPa (6.9 ± 1.4 bar, 100 ± 20 psi).

Note Air in the helium line must be purged or given sufficient time to be purged for normal performance. ▲

3. Click the **Display Status View** button in the Tune Plus window. Check whether the pressure measured by the ion gauge is $\leq 5 \times 10^{-9}$ mbar, and the pressure measured by the Pirani gauge is around 1 mbar. Compare the values of the other parameters in the status panel with values that you recorded previously.
4. In case of an Orbitrap Velos Pro ETD mass spectrometer, start up the ETD Module as described in “Starting the ETD Module After a Complete Shutdown”. In case of an Orbitrap Velos Pro mass spectrometer, continue to set up for ESI or APCI operation as described in *Orbitrap Velos Pro Getting Started*.

Starting the ETD Module After a Complete Shutdown

❖ To start up the ETD Module after a complete shutdown

1. Start the Orbitrap Velos Pro ETD mass spectrometer according to the start up procedures given in “Starting Up the System after a Shutdown” above. This also turns on the ETD Module as the ETD Module power controls are linked to the MS power controls (see “ETD Module Interface Board” on page 1-25).
2. If the Tune Plus window is not already open, choose **Start > Programs > Thermo Instruments > LTQ > LTQ Tune** from the taskbar. The Tune plus window will open.



You can determine the state of the MS detector by observing the state of the On/Off/Standby button on the Control/Scan Mode toolbar. (See Figure 2-2 on page 2-4.) The three different states of the On/Standby button are shown at the left.

3. Click the **Display Status View** button in the Tune Plus window. Check the reagent vacuum parameters:
 - Reagent ion gauge pressure: 20 to 35×10^{-5} Torr
 - Reagent Convectron gauge pressure: <0.08 Torr
 - Reagent turbomolecular pump speed : > 90%
4. Continue to set up the instrument for operation as described in *Orbitrap Velos Pro Getting Started* manual.

Resetting the System

If the communication link between Orbitrap Velos Pro mass spectrometer and data system computer is lost, it may be necessary to reset the system using the Reset button of the LTQ Velos Pro mass spectrometer.

The procedure given here assumes that the Orbitrap Velos Pro mass spectrometer and data system computer are both powered on and are operational. If the instrument, data system computer, or both are off, see [“Starting Up the System after a Shutdown”](#) on [page 2-9](#).

To reset the Orbitrap Velos Pro mass spectrometer, press the Reset button of the LTQ Velos Pro. See the *LTQ Series Hardware Manual* for the location of the Reset button. When you press the Reset button, the following occurs:

1. An interrupt on the mainboard of the internal computer causes the internal computer to reboot. All LEDs on the front panel are off except the Power LED.
2. After several seconds, the Communication LED turns yellow to indicate that the data system and the instrument are starting to establish a communication link.
3. After several more seconds, the Communication LED turns green to indicate that the data system and the instrument have established a communication link. Software for the operation of the instrument is then transferred from the data system to the instrument.
4. After three min, the software transfer is complete. The System LED turns either green to indicate that the instrument is functional and the high voltages are on or yellow to indicate that the instrument is functional and it is in Standby condition.

Resetting Tune and Calibration Parameters to their Default Values

You can reset the Orbitrap Velos Pro system tune and calibration parameters to their default values at any time. This feature may be useful if you have manually set some parameters that have resulted in less than optimum performance.

❖ To reset the Orbitrap Velos Pro MS tune and calibration parameters

In the Tune Plus window,

- Choose **File > Restore Factory Calibration** to restore the default calibration parameters, or
- Choose **File > Restore Factory Tune Method** to restore the default tune parameters.

Note Make sure that any problems you might be experiencing are not due to improper API source settings (spray voltage, sheath and auxiliary gas flow, ion transfer capillary temperature, etc.) before resetting the system parameters to their default values. ▲

Turning Off the Reagent Ion Source: What to Expect

In the Orbitrap Velos Pro ETD mass spectrometer, the reagent ion source controls can be accessed as described in “Placing the ETD Module in Standby Condition” on page 2-4. When you deselect the Reagent Ion Source On check box in the Reagent Ion Source dialog box (See Figure 2-4.), the ETD source and reagent heaters are placed in Standby condition.

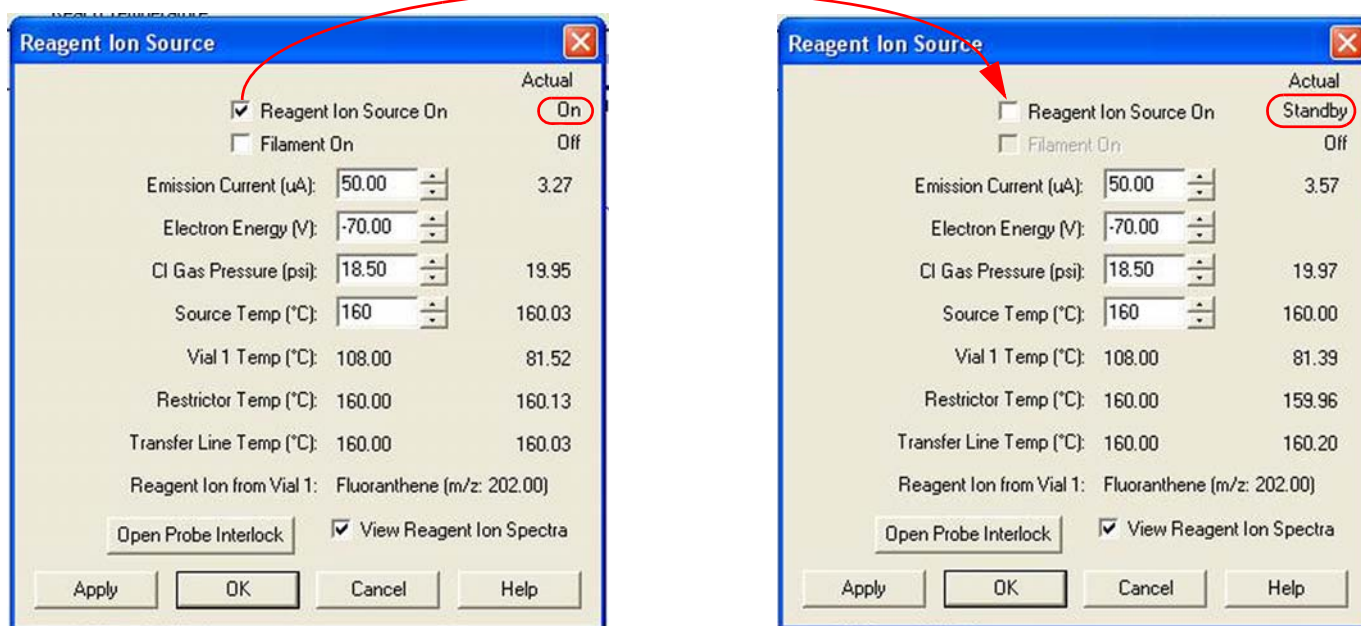


Figure 2-4. Placing the reagent ion source in Standby condition

When the ETD Module is placed in Standby condition, the filament and vial heaters are turned off. Simultaneously a valve opens to allow nitrogen gas to cool the reagent vials. This cooling nitrogen runs until the vials reach 70 °C. The audible rush (hissing noise) of nitrogen from the reagent ion source area in the back of the ETD Module is normal operation.



Warning Burn Hazard. The reagent vials are too hot to handle after the cooling nitrogen turns off at a vial temperature of 70 °C. Verify that the reagent vials have cooled down to a safe temperature before handling them. This can take up to 90 minutes after the cooling nitrogen has turned off. ▲

Other conditions that will cause the ETD Module to remain in Standby:

- Attempting to turn on the reagent ion source when the restrictor heater, transfer line heater, and the source heater are not at their target temperatures.

- Whenever either the mass spectrometer or the ETD Module goes into Standby mode. Reagent vial nitrogen cooling will turn on if the vials are at an elevated temperature.



Exception: If the Orbitrap Velos Pro ETD mass spectrometer is placed in Standby by clicking the Standby button in Tune Plus (see Standby icon in the margin), there is an hour delay before the cooling nitrogen turns on. ▲

- Whenever the pressure in the mass spectrometer or the ETD Module exceeds its protection limit. Reagent vial nitrogen cooling will turn on if the vials are at an elevated temperature.
- Whenever the abundance of reagent ions becomes insufficient as determined by the AGC setting. When this occurs, the Orbitrap Velos Pro ETD mass spectrometer completes the Xcalibur Sequence step in progress before going into Standby mode.

Chapter 3 User Maintenance

This chapter describes routine maintenance procedures that must be performed to ensure optimum performance of the Orbitrap Velos Pro mass spectrometer.

For instructions on maintaining the LTQ Velos Pro linear trap, refer to the *LTQ Series Hardware Manual*. For instructions on maintaining LCs or autosamplers, refer to the manual that comes with the LC or autosampler.

Note It is the user's responsibility to maintain the system properly by performing the system maintenance procedures on a regular basis. ▲

The following topics are described in this chapter:

- “General Remarks” on page 3-2
- “Baking Out the System” on page 3-4
- “Maintenance of the Vacuum System” on page 3-4
- “Maintenance of the ETD Module” on page 3-12
- “Maintenance of the Cooling Circuit” on page 3-58

General Remarks

Preventive maintenance must commence with installation, and must continue during the warranty period to maintain the warranty. Thermo Fisher Scientific offers maintenance and service contracts. Contact your local Thermo Fisher Scientific office for more information. Routine and infrequent maintenance procedures are listed in [Table 3-1](#).

Table 3-1. User maintenance procedures

MS Component	Procedure	Frequency	Procedure Location
Analyzer	System bakeout	If necessary (e.g. after performing maintenance work on the vacuum system)	page 3-4
Rotary-vane pumps	Add oil	If oil level is low	Manufacturer's documentation
	Change oil	Every three months or if oil is cloudy or discolored	Manufacturer's documentation page 3-5
Turbomolecular pumps	Replace operating fluid reservoir and pump bearings	Every four years	Manufacturer's documentation page 3-11
Cooling water circuit	Check cooling fluid level Check cooling fluid filter Check air inlet filter	See manufacturer's documentation	Manufacturer's documentation page 3-58
	Replace filter cartridge	Annually	page 3-58
ETD Module	Clean ion volume	As needed ^a	page 3-21
	Replace inlet valve components	As needed ^a	page 3-45
	Clean ion source lenses	As needed ^a	page 3-33
	Clean ion source	As needed ^a	page 3-40
	Replace ion source filament	As needed ^a	page 3-42
	Replace reagent vials	As needed ^a	page 3-48
	Check rotary-vane pump oil and add when needed	Every month	page 3-8
	Change rotary-vane pump oil	Every four months	page 3-9
Clean rear cooling fans	Every four months	page 3-57	

^a As needed depends on how close the component is to the electron transfer reagent introduction point. For example, the ion volume is closer to the fluoranthene introduction point than any other component and requires the most frequent cleaning.

To successfully carry out the procedures listed in this chapter, observe the following rules:

- Proceed methodically.
- Always wear clean, lint-free, and powder-free gloves when handling the components of the API source, ion optics, mass analyzer, and ion detection system.

Thermo Fisher Scientific recommends the following gloves: white nitrile clean room gloves (Fisher Scientific P/N 19-120-2947B [size medium] or P/N 19-120-2947C [size large]; Thermo Scientific P/N 23827-0008) [size medium] or P/N 23827-0009 [size large]).

- Never re-use gloves after you remove them because the surface contaminants on them will re-contaminate clean parts.
- Always wear protective eye wear when you clean parts.
- Always place the components on a clean, lint-free, and powder-free surface.
- Always cover the opening in the top of the vacuum manifold with a large, lint-free tissue whenever you remove the top cover plate of the vacuum manifold.
- Never overtighten a screw or use excessive force.
- Dirty tools can contaminate your system. Keep the tools clean and use them exclusively for maintenance and service work at the Orbitrap Velos Pro mass spectrometer.
- Never insert a test probe (for example, an oscilloscope probe) into the sockets of female cable connectors on PCBs.

Returning Parts

To protect our employees, we ask you for some special precautions when returning parts for exchange or repair to the factory. Your signature on the [Repair Covering letter](#) confirms that the returned parts have been de-contaminated and are free of hazardous materials. See ["Safety Advice for Possible Contamination"](#) on [page ix](#) for further information.

Cleaning the Surface of the Instrument

Clean the outside of the instrument with a dry cloth. For removing stains or fingerprints on the surface of the instrument (panels, for example), slightly dampen the cloth (preferably made of microfiber) with distilled water.

Caution Prevent any liquids from entering the inside of the instrument. ▲

Maintenance of the Vacuum System

This section contains instructions for performing a system bakeout and for performing pumps maintenance.

Baking Out the System

Collected or remaining gases and molecules as well as moisture can lead to an increased number of collisions with sample ions in the high vacuum region of the instrument. The bakeout procedure removes these contaminations. Therefore, we recommend to bake out the instrument if the high vacuum decreases noticeably during routine operation.

Bakeout is mandatory after maintenance or service work is performed in the analyzer region where the system is vented.

Note Pumping down the system after venting takes at least eight hours, and usually requires overnight baking of the system. ▲

In case the system has been vented during a power failure, it is necessary to bake out the system to obtain the operating vacuum. See “[Behavior of the System in Case of a Main Failure](#)” on [page 2-2](#).

Bakeout Procedure

❖ To perform a system bakeout

1. Place the system in Standby condition as described in “[Placing the Instrument in Standby Condition](#)” on [page 2-4](#).
2. Put the FT Electronics switch at the power control panel into the On position.



Figure 3-1. Bakeout timer

3. Set the bakeout time by entering the desired time (hh:mm) with the up/down keys of the bakeout timer. See [Figure 3-1](#) on [page 3-4](#).



4. Start the bakeout procedure by pressing the green start button on the right. The Orbitrap Velos Pro mass spectrometer indicates a running bakeout procedure by the flashing Vacuum and System LEDs on the front side of the instrument. See [Figure 1-4](#) on [page 1-5](#).



You can stop a running bakeout procedure by pressing the orange reset/stop button on the left side. Also press this button after you have changed the preset bakeout time.

5. The bakeout procedure is terminated because of two reasons:
 - The preset duration has expired, or
 - The vacuum has risen above a preset value.

The termination of the baking process is indicated by the status LEDs (System and Vacuum) on the front side that have stopped flashing.

Maintenance of the Forepumps

Rotary-vane pumps require minimal maintenance. All that is required to maintain the rotary-vane pump is to inspect, add, purge, and change the pump oil.

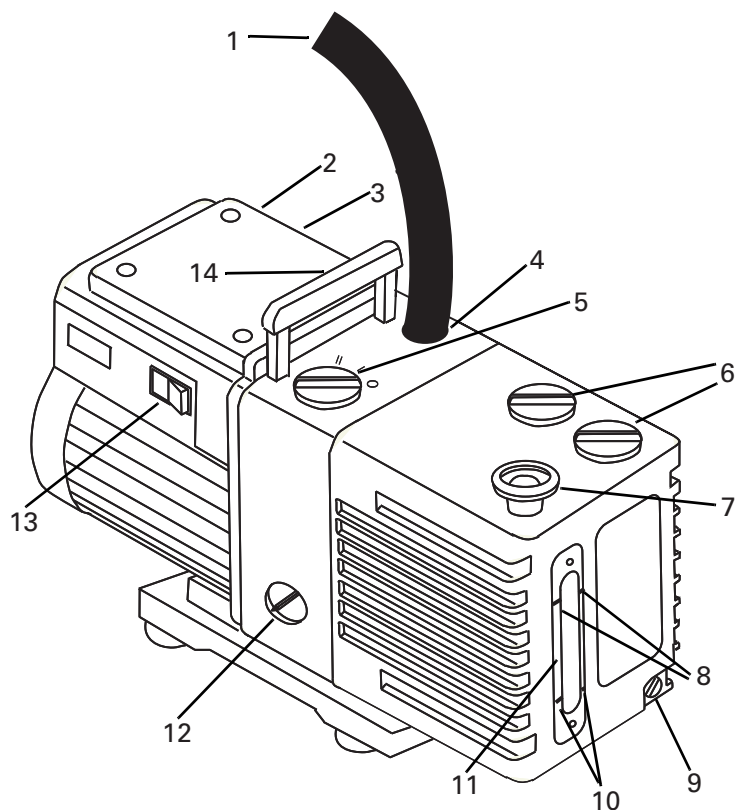
For maintenance of the forepumps of the MS portion, refer to the *LTQ Series Hardware Manual* or the pump manufacturer's manual.

Note The manuals of the pump manufacturers give detailed advice regarding safety, operation, maintenance, and installation. Please note the warnings and precautions contained in these manuals! ▲

Maintenance of the ETD Forepump

Rotary-vane pump oil (P/N A0301-15101) is a translucent light amber color and it should be checked often. During normal operation, oil must always be visible in the oil level sight glass between the MIN and MAX marks. If the oil level is below the MIN mark, add oil. If the oil is cloudy or discolored, purge the oil to decontaminate dissolved solvents. If the pump oil is still discolored, change it. You should change the pump oil every 3000 hours (about four months) of operation.

The rotary-vane pump major components are shown in [Figure 3-2](#) on [page 3-6](#).



Labeled components: 1=Foreline Vacuum Hose, 2=Electrical Inlet Connector, 3=Voltage Indicator, 4=Inlet Port, 5=Gas Ballast Control, 6=Oil Filler Plugs, 7=Outlet Port, 8=MAX Marks, 9=Oil Drain Plug, 10=MIN Marks, 11=Oil Level Sight Glass, 12=Mode Selector, 13= On/Off Switch, 14=Lifting Handle

Figure 3-2. Schematic of ETD forepump

Note During normal operation, the mode selector switch is set to high-vacuum mode (turned fully clockwise) and the gas-ballast control is closed (0). ▲

Accessing the ETD Forepump

As described in “[Forepump of the ETD Module](#)” on [page 1-34](#), the ETD forepump is located in a cabinet at the bottom of the ETD Module. To access the ETD forepump, you have to remove the lower panel as indicated in [Figure 3-3](#) on [page 3-7](#).

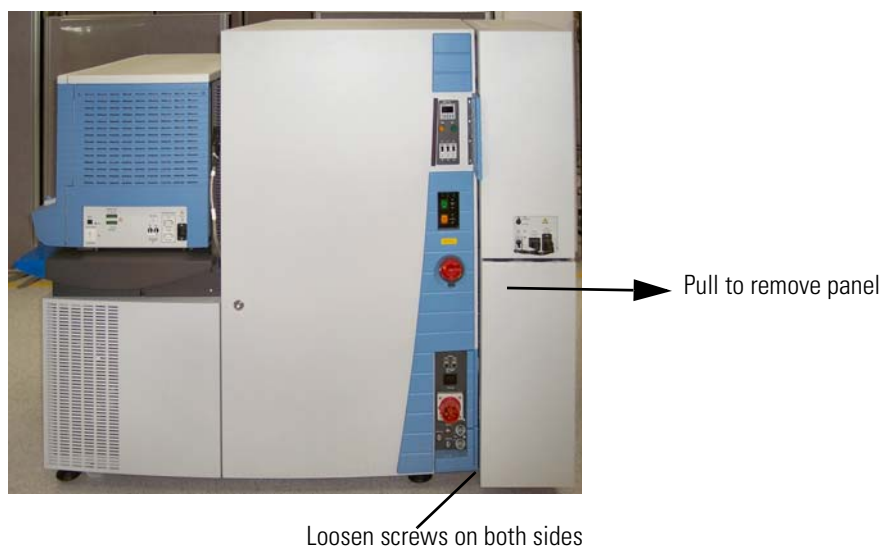


Figure 3-3. Accessing the ETD Forepump: Removing the panel

Two pairs of hooks under the top panel hold the bottom panel. They mount into corresponding openings at the top side of the bottom panel. [Figure 3-4](#) shows the details for the right side of the instrument. The panel hangs on the hooks and comes off if lifted up a little and getting pulled on into the backwards direction.

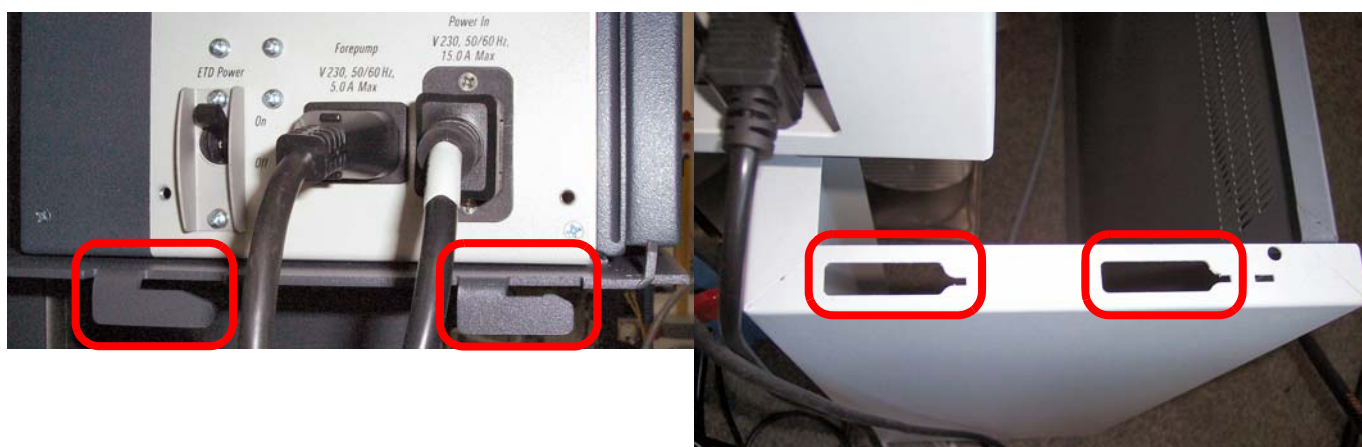


Figure 3-4. Hooks (left) and top side of detached bottom panel (right)

On the bottom of the rear side of the MS portion, two Allen screws fix the panel to the instrument frame by means of fork-like extensions (lugs). See [Figure 3-5](#) on [page 3-8](#).



Figure 3-5. Lugs for fixing the bottom panel

❖ **To remove the panel**

1. Use a 6 mm Allen wrench to loosen the screws that fix the bottom panel. Take care not to loosen the screws completely.
2. Pull the panel horizontally away from the instrument until it comes clear from the hooks.
3. Remove the panel from the instrument and store it at a safe place.

To reattach the panel, proceed in the reverse order.

Adding Oil to the ETD Forepump

The pump oil level must be between the MIN and MAX marks on the oil level sight glass for the pump to operate properly. Pump oil (P/N A0301-15101) is added as needed when the oil level is below the MIN mark on the oil level sight glass.

You can check the oil level by looking at the oil level sight glass, which is shown in [Figure 3-2](#) on [page 3-6](#). If the ETD forepump oil level is low, follow these steps to add more oil.

❖ **To add oil to the ETD forepump**

1. Shut down and vent the Orbitrap Velos Pro ETD mass spectrometer.

Caution Shut down and unplug the instrument before adding oil. ▲

2. Remove the lower panel at the rear side of the ETD Module as described on [page 3-8](#).
3. Remove one of the oil filler plugs from the rotary-vane pump.

Caution To maintain optimal performance and prevent damage to the ETD forepump, only use factory-approved rotary-vane pump oil. ▲

4. Add fresh oil to the reservoir until the oil is half way between the MIN and MAX level marks. If the oil level goes above the

MAX level mark, remove the drain plug and drain the excess oil into a suitable container.

5. Insert the oil filler plug back into the rotary-vane pump.
6. Reattach the lower panel at the rear side of the ETD Module.
7. Plug in the instrument.
8. Restart the system.

Purging the Rotary-Vane Pump Oil

When the rotary-vane pump oil becomes cloudy or discolored, purge the oil. Purging (or decontaminating) the oil removes dissolved gases and low boiling-point liquids. You can purge the oil without interrupting system operation, but do not purge it during an acquisition or while the electron multiplier or filament is powered on.

❖ To purge the rotary-vane pump oil

1. Remove the lower panel at the rear side of the ETD Module as described on [page 3-8](#).
2. Set the gas ballast control (See [Figure 3-2](#) on [page 3-6](#).) to Low Flow (I).
3. Operate the pump for 10 minutes or until the oil is clear. If the oil remains cloudy or discolored after 10 minutes, replace the oil.
4. Set the gas ballast control to Closed (0), as shown in [Figure 3-6](#).

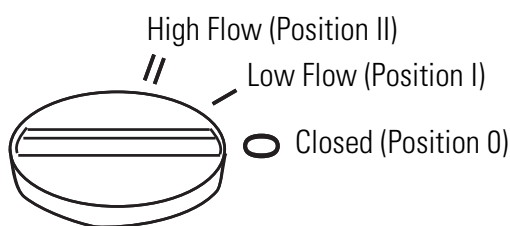


Figure 3-6. Gas ballast control positions

5. Reattach the lower panel at the rear side of the ETD Module.

Changing the Rotary-Vane Pump Oil

You should change the ETD forepump oil every four months (about 3000 hours of operation).

Supplies needed for changing the ETD forepump oil:

- Rotary-vane pump oil (P/N A0301-15101)
- Suitable container for removing spent or excess oil

Note For best results, change the oil while the ETD forepump is still warm after operation. Be careful, however, as the oil can still be very hot at this time. ▲



Warning Burn Hazard. Handle hot pump oil carefully to avoid being burned or injured. ▲

❖ **To change the ETD forepump oil**

1. Shut down and vent the Orbitrap Velos Pro ETD mass spectrometer.

Caution Shut down and unplug the instrument before adding oil. ▲

2. Remove the lower panel at the rear side of the ETD Module as described on [page 3-8](#).
3. Disassemble the rotary-vane pump.
 - a. Disconnect the foreline vacuum hose. (See [Figure 3-2](#) on [page 3-6](#).)
 - b. Disconnect the exhaust vacuum hose.
 - c. Place the rotary-vane pump on a bench.



Warning Lifting Hazard. Use the proper lifting technique to lift the ETD forepump. It weighs approximately 23 kg (50 lb). ▲

4. Drain the spent oil.
 - a. Remove one of the oil filler plugs.
 - b. Remove the oil drain plug and allow the oil to drain into a suitable container.
 - c. Dispose of the spent oil according to local environmental regulations.
 - d. Replace the oil drain plug.
5. Add fresh oil.
 - a. Add oil into oil filler reservoir half way between the MIN and MAX level marks.

- b. If the oil level goes above the MAX level mark, remove the drain plug and drain the excess oil from the pump.
6. Reassemble the rotary-vane pump.
 - a. Replace the oil filler plug.
 - b. Return the rotary-vane pump to the floor.
 - c. Reconnect the foreline vacuum hose to the rotary-vane pump.
 - d. Reconnect the exhaust vacuum hose to the rotary-vane pump.
 - e. Plug in the rotary-vane pump.
7. Reattach the lower panel at the rear side of the ETD Module.
8. Plug in the instrument.
9. Restart the system.

Maintenance of the Turbomolecular Pumps

The turbomolecular pumps in the MS portion of the Orbitrap Velos Pro mass spectrometer need maintenance work by the user that is briefly outlined below. In contrast, the turbomolecular pump in the ETD Module of the Orbitrap Velos Pro ETD mass spectrometer contains no user-serviceable parts.

Note The manuals of the pump manufacturers give detailed advice regarding safety, operation, maintenance, and installation. Please note the warnings and cautions contained in these manuals! ▲

Replacing the Operating Fluid Reservoir of the Turbomolecular Pumps

The manufacturer recommends replacing the operating fluid reservoirs of the turbomolecular pumps at least every four years. The storage stability of the operating fluid is limited. The specification of durability is given by the pump manufacturer. The disposal of used oil is subject to the relevant regulations.

Replacements for the operating fluid reservoirs including Porex rods (HiPace™ 80: P/N 1275740; HiPace™ 300: P/N 1275730) are available from Thermo Fisher Scientific.

Note The pump bearings have also to be replaced at least every four years. This maintenance operation requires special training and additional equipment. Therefore, Thermo Fisher Scientific recommends calling a Thermo Fisher Scientific field service engineer to replace both the operating fluid reservoir and the pump bearings. ▲

Maintenance of the ETD Module

This section describes routine ETD Module maintenance procedures that must be carried out to ensure optimum performance of the system. Some of the procedures describe how to clean components of the ETD Module. Others involve replacing components or replenishing the electron transfer reagent. See also the *ETD Module Hardware Manual* for additional information.

Figure 3-7 illustrates the sequence in which to perform routine maintenance on the ETD system.

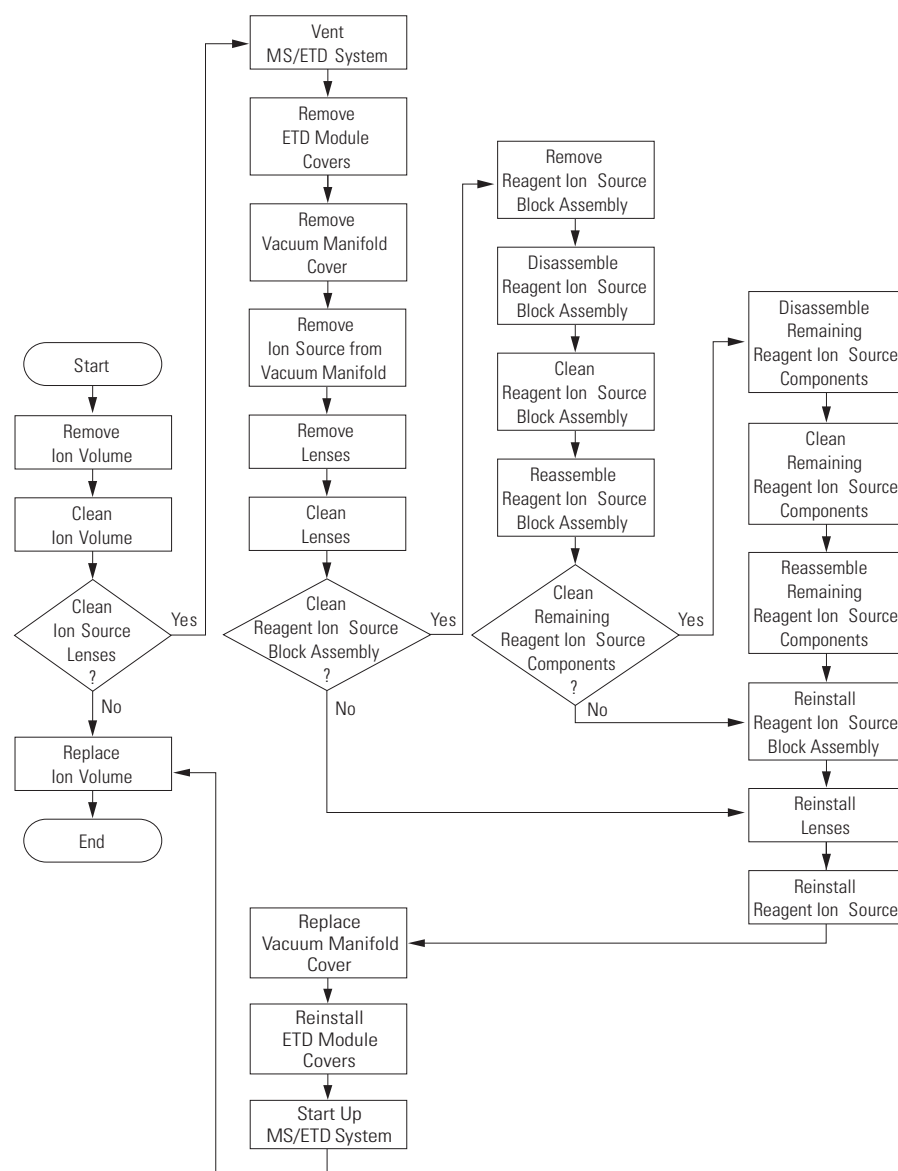


Figure 3-7. Routine maintenance sequence for ETD system

Table 3-2 on page 3-13 gives advice for correcting frequent problems with the ETD system.

Table 3-2. Indications requiring maintenance of the ETD system

Symptom	Cause	Fix
No ions at m/z 202 with the emission current at the correct level.	The m/z 202 is outside the mass range.	Set the starting mass lower.
The m/z 202 signal intensity drops slowly over several days when the emission current is at the correct level.	The ion volume needs to be cleaned or replaced.	Clean or replace the ion volume when the injection time is over 100 ms. See page 3-21 .
A system error message advising that the maximum injection time has been reached for the ETD AGC.	The AGC target has not been reached within the specified time limit. The ion volume needs to be cleaned or replaced.	Clean the ion volume. Increase the maximum injection time limit.
Sudden and complete drop of m/z 202 level, low emission current.	The filament may just have blown out.	Check the filament. Replace it if necessary. See page 3-42 .

Handling and Cleaning Reagent Ion Source Parts

A large part of maintaining your reagent ion source consists of making sure that all the components are clean. Use the cleaning procedures listed in this section to clean stainless steel and non-stainless steel parts. However, use caution when doing so, because some components can be damaged by exposure to liquids.

How often you clean the reagent ion source depends on the amount of reagent introduced into the system. In general, the closer a component is to where the reagent ion is introduced, the more rapidly it becomes dirty (see the footnote in [Table 3-1](#) on [page 3-2](#)). For example, the ion volume needs to be cleaned more often than other parts.

Many parts can be removed and disassembled by hand. Make sure you have all the necessary tools before carrying out a procedure. See below for a list of the tools and supplies generally needed for maintenance of the reagent ion source. Tools should be used only for the maintenance of the reagent ion source and be free of grease or other residues. Handle parts in a manner that maintains their cleanliness.

Note It is crucial that the cleanliness of the parts be maintained when they are handled. Wear gloves and place the parts on surfaces that are clean if the parts are not returned directly to the instrument. If clean surfaces are not available, place the parts on fresh lint free wipes or cloths or aluminum foil that has not been used for any other purpose. ▲

The following tools and supplies are needed for reagent ion source maintenance:

- Clean, dry gas (air or nitrogen)
- New, white nitrile clean room gloves (Fisher Scientific P/N 19-120-2947B [size medium] or P/N 19-120-2947C [size large]; Thermo Scientific P/N 23827-0008) [size medium] or P/N 23827-0009 [size large])
- Lint-free cloth or paper
- Nut driver, 5.5 mm
- Protective eyewear
- Screwdriver, Phillips #2
- Screwdriver, flat blade
- Wrench, adjustable
- Wrench, Allen, 2 mm, 2.5 mm, 3 mm, 4 mm, 5/32-in., 5/64-in., 1/16-in
- Wrench, open-ended, 1/4-in., 5/16-in., 7/16-in. (2), 1/2-in., 9/16-in.
- Wrench, socket, 1/2-in.

Cleaning Stainless Steel Parts

The reagent ion source, ion volume assembly, ion source block, and lenses are made from stainless steel. To clean these parts, follow the procedure described in this topic. Use this procedure with caution because some components can be damaged when exposed to liquids.

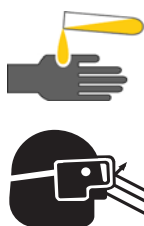
The following tools and supplies are needed for cleaning stainless steel parts in the reagent ion source:

- Acetone, analytical grade (or other suitable solvent)
- Aluminum oxide abrasive, number 600 (P/N 32000-60340)
- Applicators, cotton-tipped
- Beaker, 450 mL
- Clean, dry gas
- De-ionized water
- Detergent (Alconox®, Micro, or equivalent)
- Dremel® rotary tool or equivalent (recommended)

- Foil, aluminum
- Forceps
- New, white nitrile clean room gloves
- Glycerol, reagent grade
- Lint-free cloth
- Protective eyewear
- Tap water
- Toothbrush, soft
- Ultrasonic cleaner

Caution Do not use this procedure to clean ceramic, aluminum, or gold plated parts. See page 3-17 for the method for cleaning ceramic, aluminum, or gold plated parts. ▲

Caution Follow the subsequent instructions precisely. If done wrong, the cleaning procedure could damage the ion source lenses. ▲



Warning Hand and Eye Hazard. Wear impermeable laboratory gloves and eye protection when performing these cleaning procedures. ▲

❖ **To clean reagent ion source stainless steel parts**

1. Remove contamination from the surfaces being cleaned.
 - a. Use a slurry of number 600 aluminum oxide in glycerol and a cleaning brush or cotton-tipped applicator. Contamination often appears as dark or discolored areas, but may not be visible. The heaviest contamination is usually found around the apertures, such as the electron entrance hole on the ion volume.
 - b. Clean each part thoroughly, even if no contamination is visible.
 - c. Use the wooden end of an applicator that is cut at an angle to clean the inside corners.
 - d. Use a Dremel® tool with the polishing swab at its lowest speed. This will increase the cleaning efficiency and decrease the time required to clean the part.



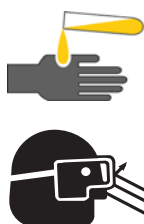
Warning Injury Hazard. To prevent personal injury, be sure to keep the Dremel tool away from possible hazards, such as standing water or flammable solvents. ▲

2. Rinse the parts with clean water.
3. Use a clean applicator or toothbrush to remove the aluminum oxide slurry. Do not let the slurry dry on the metal because dried aluminum oxide is difficult to remove.
4. Place the parts in a warm detergent solution in an ultrasonic bath and sonicate them.
 - a. Make a solution of detergent and warm tap water in a 400 mL glass beaker.
 - b. Using forceps, place the parts in a beaker containing the warm detergent solution.
 - c. Place the beaker and contents in an ultrasonic bath for five minutes.
 - d. Rinse the parts with tap water to remove the detergent.
5. Sonicate the parts in deionized water.
 - a. Using forceps, place the parts in a beaker containing deionized water.
 - b. Place the beaker and contents in an ultrasonic bath for five minutes.
 - c. If the water is cloudy after sonicating, pour off the water, add fresh water, and place the beaker and its contents in a ultrasonic bath again for five minutes. Repeat until the water is clear.
6. Sonicate the parts in acetone.
 - a. Using forceps, place the parts in a beaker containing fresh acetone.
 - b. Place the beaker and contents in an ultrasonic bath again for five minutes.
7. Blow-dry the parts immediately. Use clean, dry gas (air or nitrogen) to blow the acetone off the parts.
8. Complete the drying process, doing one of the following:
 - Using forceps, place the parts in a 500 mL glass beaker, cover the beaker with aluminum foil, and put the beaker in an oven set at 100 °C for 30 minutes.

- Lay the parts on clean aluminum foil (dull side up) and allow to dry for 30 minutes.
9. Allow the parts to cool before reassembling them.

Cleaning Non-Stainless Steel or Hybrid Parts

To clean the stainless-steel portion of hybrid parts, follow [step 1](#) and [step 2](#) of the instructions on [page 3-15](#). Perform these steps only on the stainless-steel surfaces of hybrid parts. Do not allow the aluminum oxide slurry to contact the aluminum, ceramic, or gold plated portions of these parts.



Warning Hand and Eye Hazard. Wear impermeable laboratory gloves and eye protection when performing these cleaning procedures. ▲

The reagent ion source heater ring, filament spacer, lens holder, and spacers are non-stainless steel parts that are made from aluminum, ceramic, or are gold plated.

❖ To clean the non-stainless-steel portions of hybrid parts

1. Scrub all of the parts with a warm detergent solution.
 - a. Make a solution of detergent and warm tap water in a 500 mL glass beaker.
 - b. Dip a clean cotton-tipped applicator in the detergent mixture and use the applicator to scrub the parts.

Note Do not soak or sonicate the parts in detergent. ▲

- c. Using forceps, rinse the parts thoroughly with tap water to remove the detergent.

Caution Do not leave aluminum parts, such as the heater ring, in the detergent. Basic solutions, like detergent, damage the surface of aluminum. ▲

2. Rinse the parts in deionized water. Using forceps, dip the parts in a beaker of deionized water. Change the water if it becomes cloudy. Do not soak or sonicate the parts.

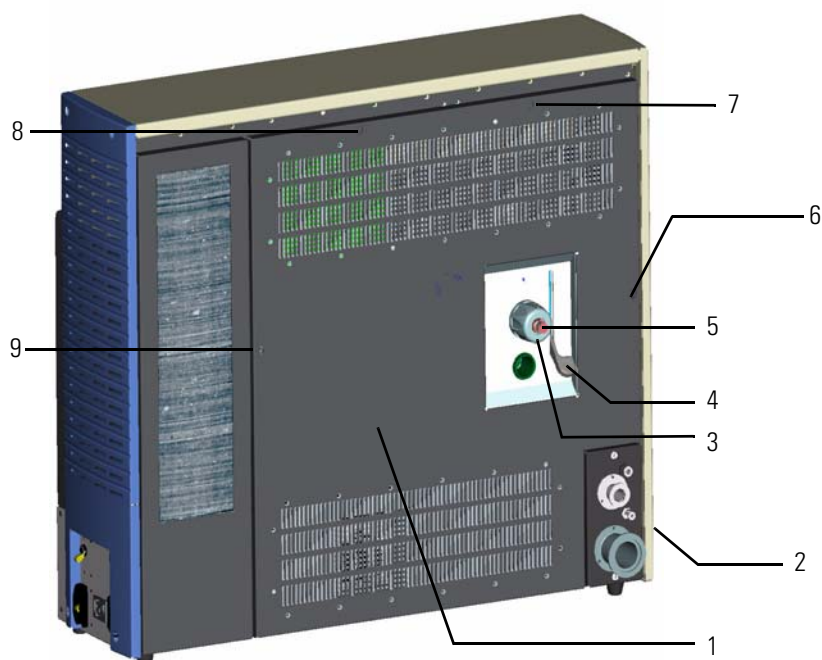
User Maintenance

Maintenance of the ETD Module

3. Rinse the parts with acetone. Using forceps, dip the parts in a beaker of acetone. Change the acetone if it becomes cloudy. Do not soak or sonicate the parts.
4. Blow-dry the parts immediately. Use clean, dry gas (air or nitrogen) to blow the acetone off the parts.

Removing the Access Panels

During some ETD Module maintenance activities, it is necessary to remove either the ETD main access panel, or the side access panel, or both (see [Figure 3-8](#)). Follow the subsequent procedures to remove these panels.



Labeled components: 1=ETD main access panel, 2=side access panel, 3=inlet valve knob, 4=inlet valve lever (down is closed, up is open), 5=inlet valve plug, 6, 7, 8, 9=panel fasteners.

Figure 3-8. Rear view of the ETD Module

Removing the ETD Main Access Panel

❖ To remove the ETD main access panel

1. Place the ETD Module to Service mode as directed in [“Placing the Instrument in Off Condition and Service Mode”](#) on [page 3-48](#).

Note In Service mode, all power to the Orbitrap Velos Pro ETD MS electronics is turned off. There are no user accessible components that carry a voltage in this mode. However, the vacuum pumps continue to operate. ▲



Warning Burn Hazard. The reagent vial heaters can be 108 °C (or set point); the transfer line, the restrictor, and the ion source can be at 160 °C. These components may be too hot to touch. Verify that all of these components are safe to touch before handling them. ▲

Note The ETD main access panel is interlocked with the ETD Module power. When the ETD main access panel is removed, all power to the ETD Module will be turned off. However, the mechanical pump and turbo pump will continue operating. ▲

2. Remove the inlet valve lever (item 4 in [Figure 3-8](#) on [page 3-18](#)) by pulling it down and away from the ETD Module main access panel. Do not rotate the lever upwards. It must remain in its down (closed) position to avoid catastrophic venting of the system.

Caution Rotating the inlet valve lever upwards (to the open position) without the inlet valve plug (item 5 in [Figure 3-8](#)) or the ion volume tool in place will cause a catastrophic venting of the system. ▲

3. Unscrew the inlet valve knob (item 3 in [Figure 3-8](#)) and remove the inlet valve plug (item 5 in [Figure 3-8](#)), the inlet valve knob, and the internal ferrule.
4. Loosen the four panel fasteners (items 6, 7, 8, and 9 in [Figure 3-8](#)).
5. The top panel rests on hooks pointing into the upward direction. Tilt the top panel towards you and lift it up and away from the ETD Module.

Removing the ETD Side Access Panel

❖ To remove the ETD side access panel

1. If it is not already in Service mode, place the ETD Module in Service mode as directed in [“Placing the Instrument in Off Condition and Service Mode”](#) on [page 3-48](#).

Note In Service mode, all power to the Orbitrap Velos Pro ETD MS electronics is turned off. There are no user accessible components that carry a voltage in this mode. However, the vacuum pumps continue to operate. ▲



Warning Burn Hazard. The reagent vial heaters can be at 108 °C (or set point); the flow restrictor, the transfer line heaters, and the ion source heater can be at 160 °C. These components may be too hot to touch. Verify that all of these components are safe to touch before handling them. ▲

Note The ETD side access panel is interlocked with the ETD Module power. When the ETD side access panel is removed, all power to the ETD Module will be turned off. However, mechanical pump and turbomolecular pump will continue operating. ▲

2. Using an Allen wrench, loosen the captive screws at the top and remove the screws at the bottom of gray plastic side panel and remove the panel.
3. Using a #2 Phillips screwdriver, loosen the three captive screws on the metal side access panel (item 2 in [Figure 3-8](#) on [page 3-18](#)) and remove the panel.



Warning Burn Hazard. Reagent vial heaters, ion source heater, flow restrictor, and transfer lines are accessible under the ETD side access panel. These are heated components. Verify that they are safe to touch before handling them. ▲

Replace the panels by following the above steps in reverse order and reversing the instructions in each step.

Maintenance of the Reagent Ion Source

The reagent ion source consists of an ion volume, filament, and ion source lenses. Because the ion volume is exposed directly to samples introduced into the reagent ion source, it requires the most frequent cleaning. You can access the ion volume assembly with or without an inlet valve.

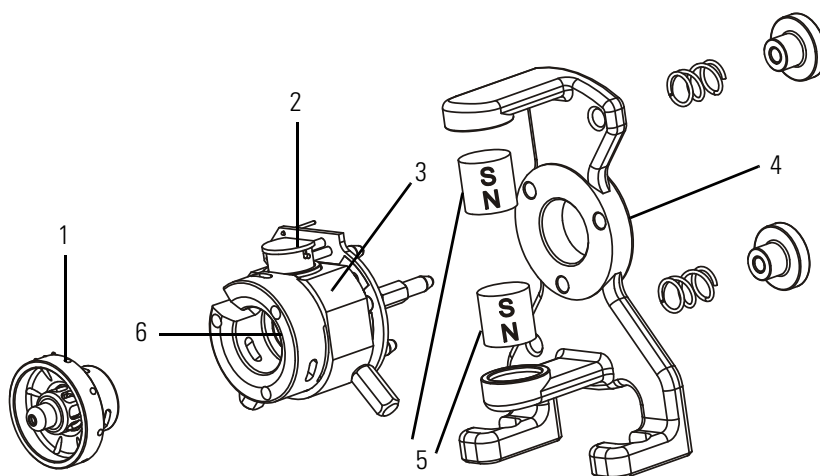
To restore system performance, always clean the ion volume first, then the ion source lenses. If cleaning either of these components does not restore system performance, try cleaning the entire reagent ion source.

This section contains these maintenance procedures:

- “Cleaning the Ion Volume With an Inlet Valve” on [page 3-21](#)
- “Cleaning the Ion Source Lens Assembly” on [page 3-33](#)
- “Cleaning the Ion Source Block” on [page 3-40](#)
- “Replacing the Ion Source Filament” on [page 3-42](#)

- “Replacing Inlet Valve Components” on page 3-45

The ion source, the ion trap, and their components are shown in Figure 3-9.



Labeled components: 1=ion source lenses, 2=filament assembly, 3=ion source block, 4=magnet support, 5=magnets, 6=ion volume (inside the ion source block, 3)

Figure 3-9. Ion source components (left view)

Cleaning the Ion Volume With an Inlet Valve

The ion volume is where molecules interact with energetic electrons to form ions. Because the ion volume is exposed directly to reagents introduced into the reagent ion source, you will have to clean it more frequently than other parts. How often you have to clean the ion volume assembly will depend on the types and amounts of reagents used.

For cleaning the ion volume with an inlet valve, the following tools and supplies are needed:

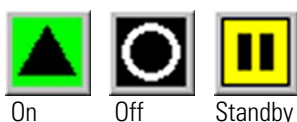
- Cleaning supplies for stainless steel parts (See “Cleaning Stainless Steel Parts” on page 3-14.)
- Gloves (clean, lint-free, and powder-free)
- Ion volume tool and guide bar
- Lint-free cloth

Using the ion volume tool allows you to access the ion volume by entering the vacuum manifold through the inlet valve without venting the instrument.

User Maintenance

Maintenance of the ETD Module

❖ To clean the ion volume with an inlet valve



1. Click the On/Standby button in the Tune Plus window to place the Orbitrap Velos Pro ETD mass spectrometer in Standby mode. See [Figure 3-10](#).

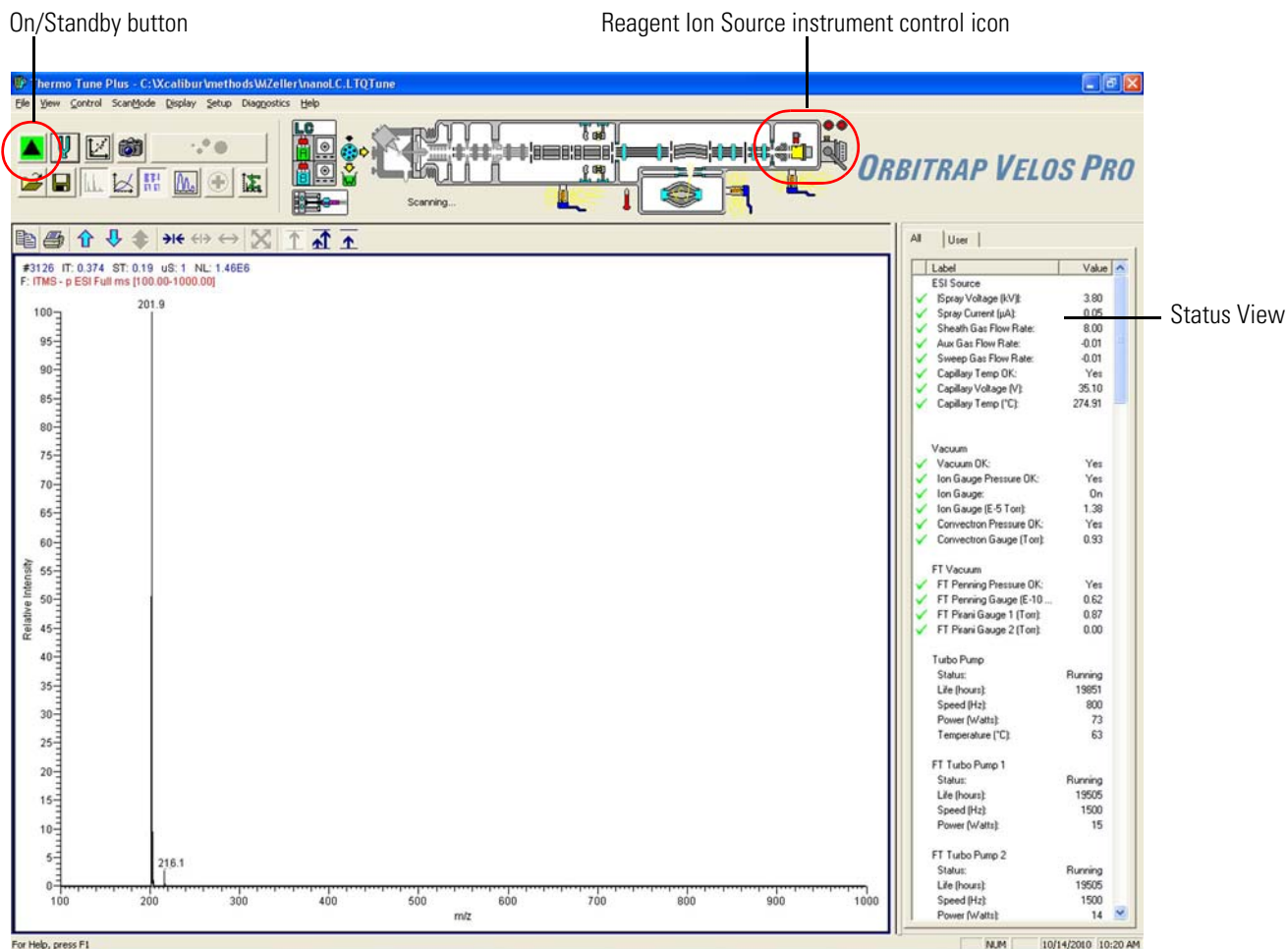
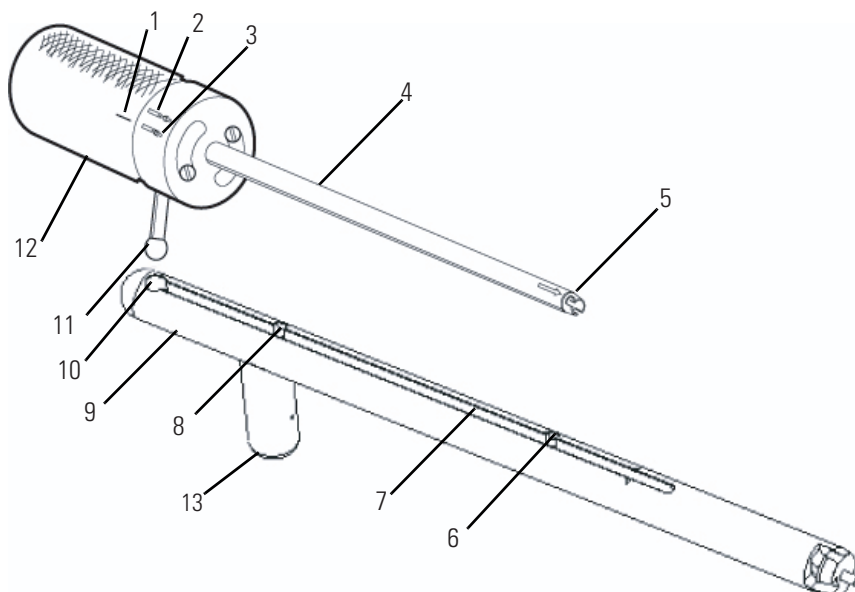


Figure 3-10. Tune Plus window (Orbitrap Velos Pro ETD)



2. Open the Reagent Ion Source dialog box ([Figure 3-17 on page 3-26](#)) in Tune Plus by clicking the Reagent Ion Source instrument control icon.
3. Place the guide bar handle (item 13 in [Figure 3-11 on page 3-23](#)) to the 3 o'clock position ([Figure 3-12 on page 3-23](#)).



Labeled components: 1=alignment line, 2=lock position, 3=unlock position, 4=ion volume tool, 5=bayonet lock, 6=second stop, 7=guide ball track, 8=first stop, 9=guide bar, 10=guide ball hole, 11=guide ball, 12=ion volume tool handle, 13=guide bar handle

Figure 3-11. Ion volume tool components

4. Insert the guide bar (item 9 in [Figure 3-11](#)) into the guide bar opening in the back of the ETD Module ([Figure 3-12](#)).

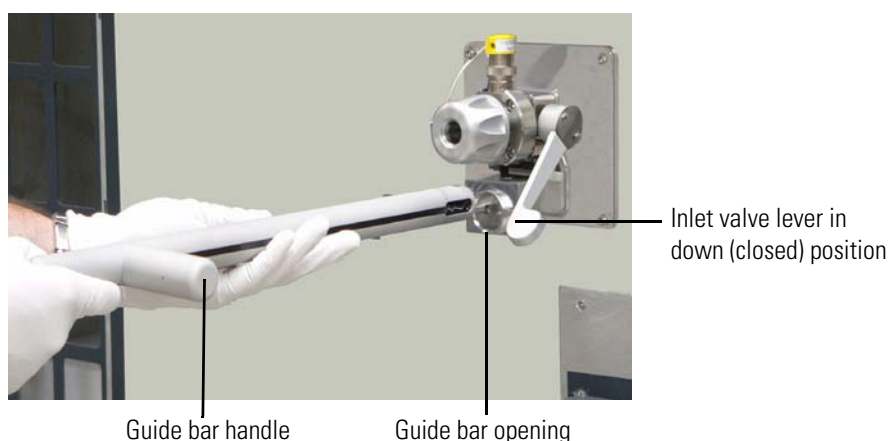


Figure 3-12. Guide bar being inserted into guide bar opening^a

^aGuide bar handle is facing to the right. The inlet valve is closed when the inlet valve lever is in the down position and open when it is in the up position.

5. Push the guide bar in as far as it will go, then rotate it 90° clockwise to lock in the guide bar ([Figure 3-13](#) on [page 3-24](#)). The guide bar handle faces the floor at the completion of this step.

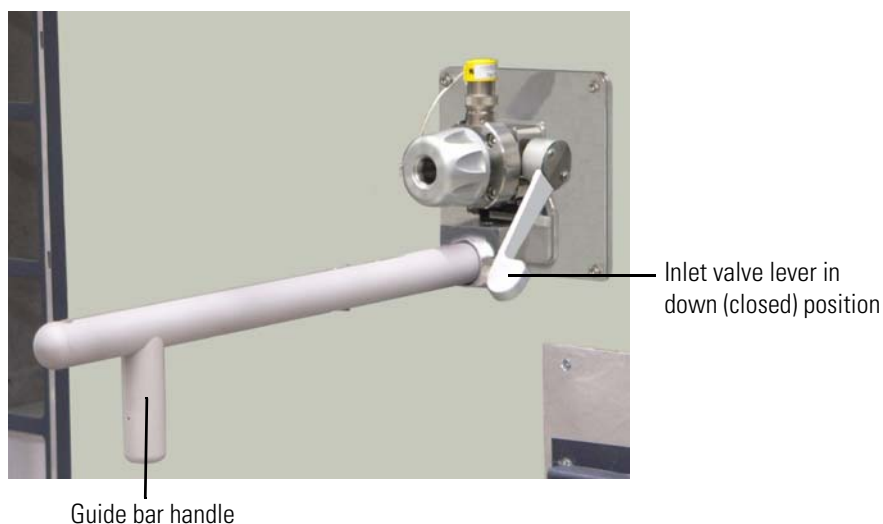
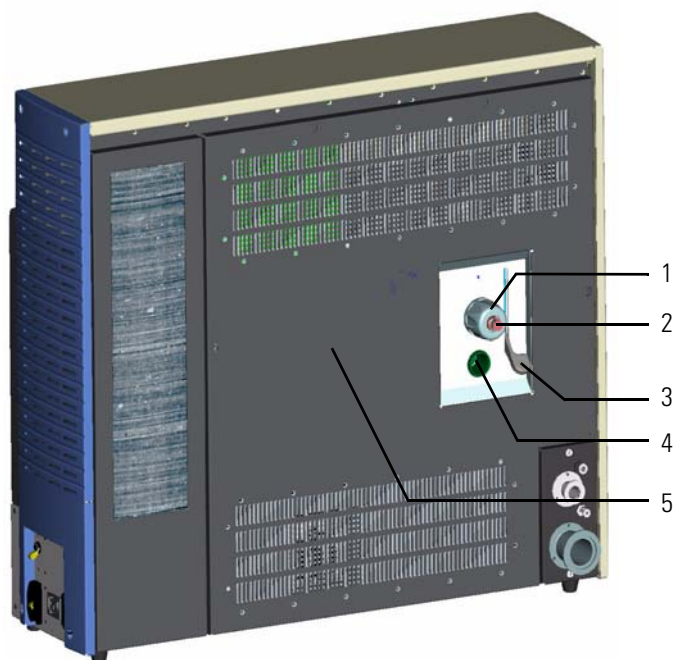


Figure 3-13. Guide bar insertion complete^a

^aGuide bar handle is facing the floor. The inlet valve is closed when the inlet valve lever is in the down position and open when it is in the up position.

6. Prepare the inlet valve and ion volume tool for insertion.



Labeled components: 1=inlet valve knob, 2=inlet valve plug, 3=inlet valve lever (down is closed, up is open), 4= guide bar opening, 5=main access panel

Figure 3-14. Rear view of the ETD Module, showing the inlet valve

Make sure the inlet valve is closed (inlet valve lever is down, as shown in [Figure 3-13](#)) and remove the inlet valve plug (item 2 in [Figure 3-14](#)). Do this by rotating (loosening) the inlet valve knob

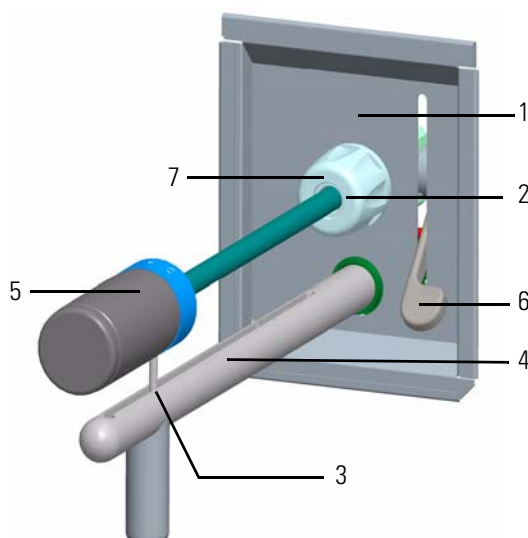
(item 1 in [Figure 3-14](#) on [page 3-24](#)) until the inlet valve plug will slide out easily. The inlet valve plug prevents air from entering the vacuum manifold in case the inlet valve is inadvertently opened.

7. Turn the ion volume tool handle to the unlock position, which indicates that the ion volume tool is in position to accept the ion volume. See [Figure 3-15](#).



Figure 3-15. Ion volume tool handle in the unlock position

8. Insert the ion volume tool and evacuate the inlet valve:
 - a. Insert the guide ball into the guide ball hole.



Labeled components: 1=ion volume tool entry housing, 2=inlet valve opening, 3=first stop, 4=guide bar, 5=ion volume tool, 6=inlet valve lever, 7=inlet valve knob

Figure 3-16. Ion volume tool guide bar first stop

- b. Slide the ion volume tool forward in the guide bar track until the guide ball is at the guide bar's first stop, which is shown in [Figure 3-11](#) on [page 3-23](#) and [Figure 3-16](#) on [page 3-25](#).
- c. Slide the ion volume tool so the guide ball is in the groove at the first stop ([Figure 3-11](#) and [Figure 3-16](#)). This prevents the probe from being pulled forward when the inlet valve is evacuated.
- d. Tighten the inlet valve knob ([Figure 3-16](#)) to ensure that a leak-tight seal is made.
- e. Click **Open Probe Interlock** in the Reagent Ion Source dialog box ([Figure 3-17](#)). A message box appears stating that the probe interlock is being pumped down. The target pressure is <0.1 mTorr. If a pressure of 0.1 mTorr or less is not obtained, replace the inlet valve seal as described in [“Replacing Inlet Valve Components”](#) on [page 3-45](#). When the target pressure is achieved, a message appears stating that the ball valve can be opened ([Figure 3-18](#)).

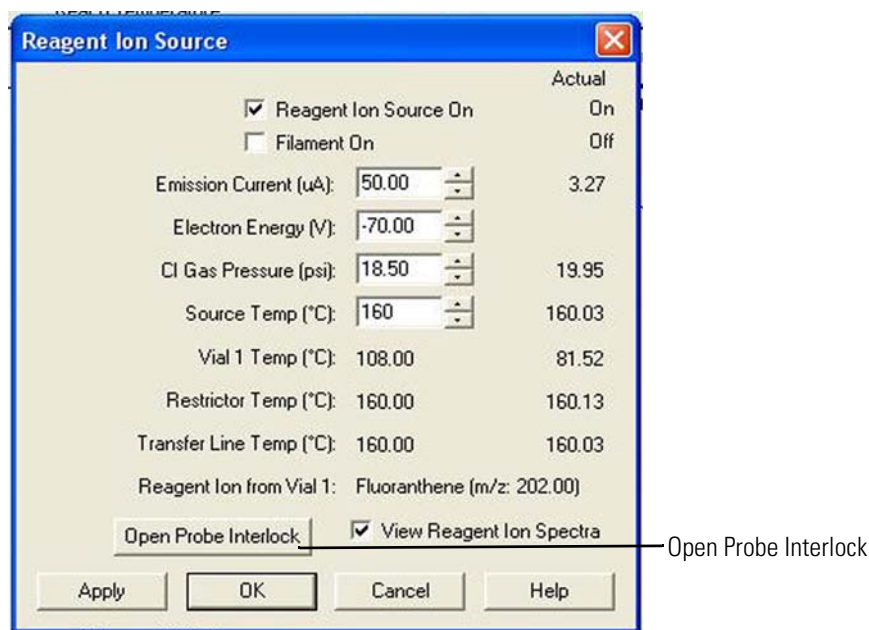
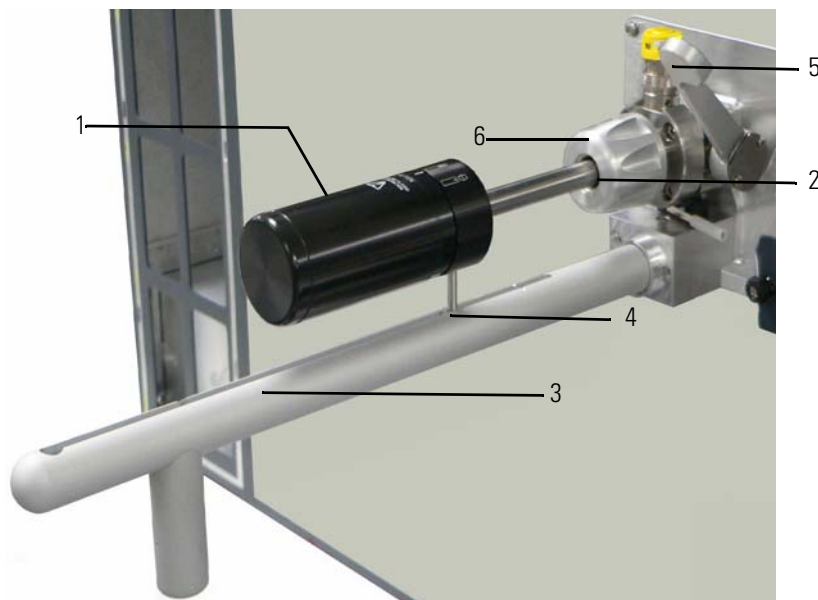


Figure 3-17. Reagent Ion Source dialog box, Open Probe Interlock button



Figure 3-18. Instrument Message box: The Ball Valve can now be opened

- f. Once evacuation is complete, push up the inlet valve lever to open the inlet valve (Figure 3-19).
9. Remove the ion volume:
- a. Slide the ion volume tool into the vacuum manifold until the tip of the ion volume tool is fully inserted into the ion volume holder, as shown in Figure 3-19.



Labeled components: 1=ion volume tool, 2=inlet valve opening, 3=guide bar, 4=second stop, 5=inlet valve lever in open (up) position, 6=inlet valve knob

Figure 3-19. Ion volume tool inserted into the inlet valve

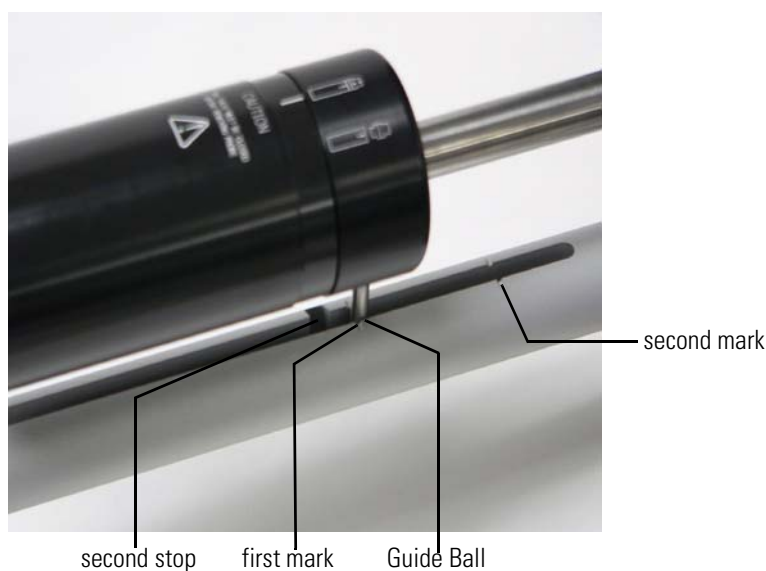


Figure 3-20. Detail of ion volume tool fully inserted into the inlet valve

You will know that the ion volume tool is fully inserted into the ion volume holder because the guide ball (item #11, [Figure 3-11](#) on [page 3-23](#)) will be just past the first mark on the guide bar as shown in [Figure 3-20](#) on [page 3-27](#).

- b. Turn the ion volume tool handle counterclockwise to the lock position, See [Figure 3-21](#). Listen for a click indicating that the handle is fully engaged in the lock position and is holding the ion volume.



Figure 3-21. Ion volume tool handle in the locked position

- c. Withdraw the ion volume tool (the ion volume is attached) until the guide ball reaches the first stop (see [Figure 3-11](#) on [page 3-23](#) and [Figure 3-16](#) on [page 3-25](#) for the first stop position).
- d. Close the inlet valve by pushing the lever down.

Caution Do not withdraw the ion volume tool beyond the point where the guide ball reaches the first stop in the guide bar. Close the inlet valve before withdrawing the ion volume tool past the first stop. Otherwise, the system will vent to the atmosphere and cleaning the components that are under vacuum will be required. ▲

- e. Loosen the inlet valve knob ([Figure 3-19](#) on [page 3-27](#)).
- f. Continue withdrawing the ion volume tool completely from the inlet valve by sliding the ion volume tool through the guide ball track in the guide bar.



Warning Burn Hazard. The ion volume will be too hot to touch. Let it cool to room temperature before handling it. ▲

10. Clean the ion volume:



- a. Turn the ion volume tool handle to the unlock position (Figure 3-15 on page 3-25). The ion volume tool handle unlock position icon is shown at the left.
- b. Remove the ion volume from the ion volume tool. Using clean gloves, press the ion volume into the tip of the ion volume tool and rotate it to disconnect the bayonet pins from the pin guides. Pull the ion volume out of the ion volume tool, as illustrated in Figure 3-22.

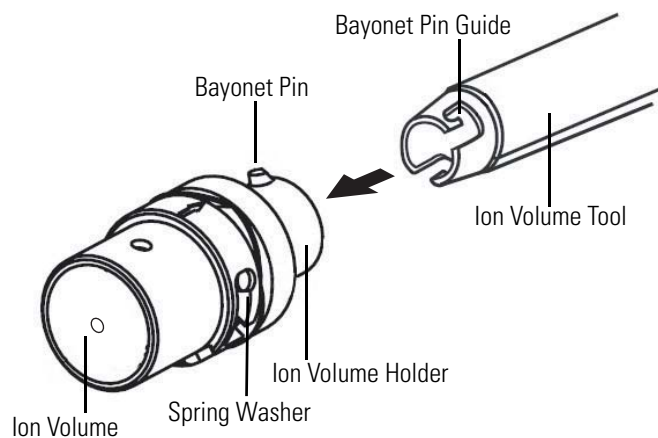


Figure 3-22. Ion volume assembly

- c. Press the ion volume into the ion volume holder and rotate the ion volume to remove it from the ion volume holder See Figure 3-23.

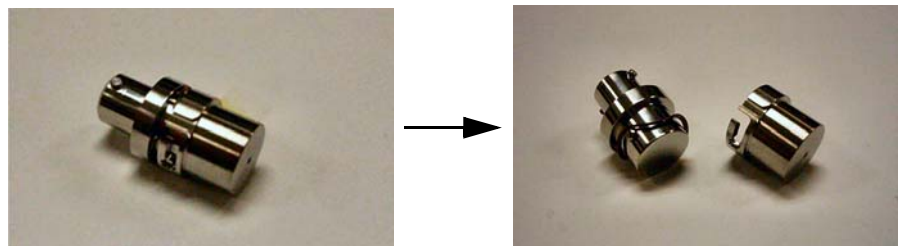


Figure 3-23. Separating ion volume and ion volume holder

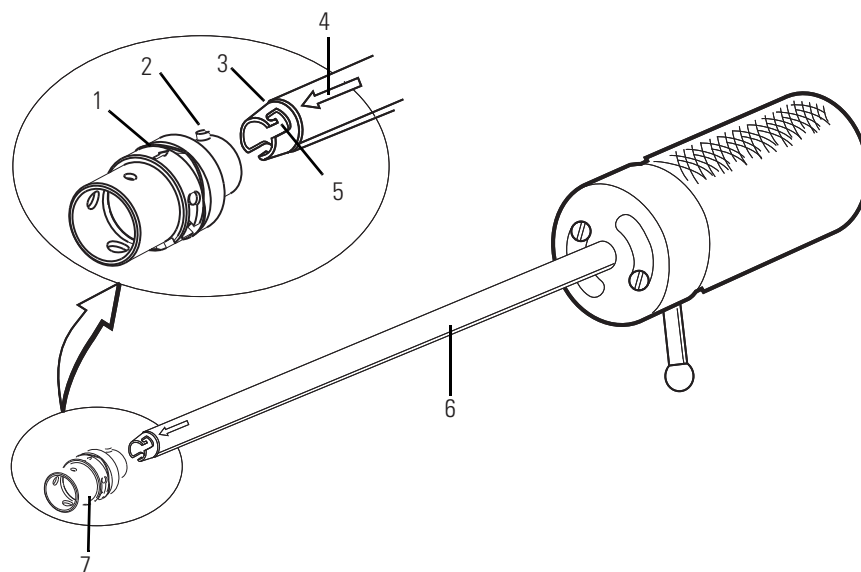
- d. Clean ion volume and ion volume holder according to the instructions in “Cleaning Stainless Steel Parts” on page 3-14.

❖ **To reinsert the ion volume**

1. Press the ion volume into the ion volume holder and rotate the ion volume to secure it to the ion volume holder.

2. Place the clean ion volume on the ion volume tool:
 - a. Place the ion volume into the bayonet lock located on the ion volume tool. Make sure that the alignment arrows on the ion volume and ion volume tool are facing each other. See [Figure 3-24](#).

Caution To avoid damage to the ion source, ensure that the arrows on the ion volume tool and ion volume are aligned. ▲



Labeled components: 1=ion volume alignment arrow, 2=bayonet pin, 3=bayonet lock, 4=ion volume tool alignment arrow, 5=bayonet guide, 6=ion volume tool, 7=ion volume

Figure 3-24. Placing the ion volume on the ion volume tool

Note Wear clean, lint-free, and powder-free gloves when you handle parts inside the vacuum manifold. ▲

- b. Turn the ion volume tool handle to the lock position. (See [Figure 3-21](#) on [page 3-28](#).)
3. Insert the ion volume tool and evacuate the inlet valve:
 - a. Insert the guide ball into the guide ball hole and slide the ion volume tool forward in the guide bar track until the guide ball is at the guide bar's first stop (see [Figure 3-11](#) on [page 3-23](#) and [Figure 3-16](#) on [page 3-25](#)).
 - b. Turn the ion volume tool so that the guide ball is in the groove at the first stop ([Figure 3-16](#) on [page 3-25](#)). This prevents the

probe from being pulled forward when the inlet valve is evacuated.

- c. Tighten the inlet valve knob to ensure a leak-tight seal (Figure 3-19 on page 3-27).
 - d. Click **Open Probe Interlock** in the Reagent Ion Source dialog box (Figure 3-17 on page 3-26). A message box will appear stating that the probe interlock is being pumped down. The target pressure is <0.1 mTorr. If a pressure of 0.1 mTorr or less is not obtained, the inlet valve seal must be replaced as described in “Replacing Inlet Valve Components” on page 3-45. When the target pressure is achieved, a message will appear stating that the ball valve can be opened. See Figure 3-18 on page 3-26.
 - e. Once evacuation is complete, push the inlet valve lever up to open the inlet valve. See Figure 3-19 on page 3-27.
4. Reinsert the ion volume:
- a. Slide the ion volume tool into the vacuum manifold, as illustrated in Figure 3-19.
 - b. Listen for a click indicating that the ion volume has connected with the ion source block. The guide ball will be slightly beyond the second stop on the guide bar. See Figure 3-20 on page 3-27.
 - c. Turn the ion volume tool handle to the unlock position. See Figure 3-25.



Figure 3-25. Ion volume tool handle in the unlock position

- i. Withdraw the ion volume tool away from the ion volume about 2.5 cm (1 in) and turn the ion volume tool handle to the lock position. See Figure 3-26 on page 3-32.
- ii. Slide the ion volume tool back into the vacuum manifold until the end of the ion volume tool just touches the ion volume.

- iii. If the ion volume tool does not go into the inlet valve completely, the ion volume is not seated properly.
- d. Withdraw the ion volume tool until the guide ball reaches the first stop (see [Figure 3-11](#) on [page 3-23](#) and [Figure 3-16](#) on [page 3-25](#)).



Figure 3-26. Ion volume tool handle in the locked position

- e. Close the inlet valve by pushing down on the inlet valve lever ([Figure 3-14](#) on [page 3-24](#)).

Caution Do not withdraw the ion volume tool beyond the point where the guide ball reaches the first stop in the guide bar. Close the inlet valve before withdrawing the ion volume tool past the first stop. Otherwise, the system vents to the atmosphere. ▲

- f. Loosen the inlet valve knob (item 6 in [Figure 3-19](#) on [page 3-27](#)).
 - g. Continue withdrawing the ion volume tool completely from the inlet valve by sliding the ion volume tool through the guide ball track in the guide bar.
5. Remove the ion volume tool and guide bar from the vacuum manifold:
 - a. Remove the guide bar by rotating it 90 degrees counter-clockwise and sliding it out of the entry housing.
 - b. Replace the inlet valve plug and tighten the inlet valve knob (item 6 in [Figure 3-19](#) on [page 3-27](#)).
 - c. Click **Close** in the message stating that the ball valve can be opened. (See [Figure 3-18](#) on [page 3-26](#).)
 6. Re-tune the MS detector.

Note Tune Plus provides an evaluation procedure for CI gas pressure under **Diagnostics > Diagnostics > Tools > System evaluation > Reagent CI gas pressure evaluation**. Thermo Fisher Scientific recommends performing this procedure after replacing the filament and/or the ion volume. ▲

Cleaning the Ion Source Lens Assembly

If cleaning the ion volume did not restore system performance, try cleaning the ion source lens assembly. The ion source lens assembly comes in direct contact with reagent ions introduced into the ETD Module and needs to be cleaned periodically (though not as often as the ion volume).

❖ To clean the ion source lens assembly

1. Prepare the ETD Module for maintenance:
 - a. Prepare a clean work area by covering the area with a clean lint-free cloth.
 - b. Shut down and vent the ETD Module. (See “[Shutting Down the Instrument](#)” on [page 2-7](#).)

Caution Shut down and unplug the Orbitrap Velos Pro ETD mass spectrometer before proceeding with the next steps of this procedure. ▲

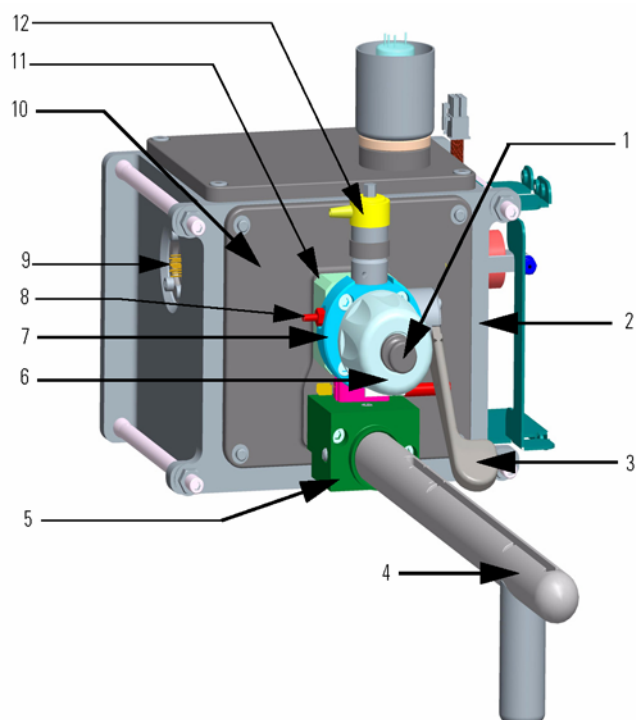
Note Wear clean, lint- and powder-free gloves when you handle parts inside the vacuum manifold. ▲



Warning Burn Hazard. The ion source may be too hot to touch even if the cooling nitrogen has completed its cycle. Be sure that the ion source has cooled to room temperature before handling it. ▲

2. Remove the ion source assembly:
 - a. Remove the main access panel of the ETD Module (item 1 in [Figure 3-8](#) on [page 3-18](#)). Follow the procedures in “[Removing the ETD Main Access Panel](#)” on [page 3-18](#).

Caution It is good practice to keep the inlet valve lever in the down (closed) position whenever it is not explicitly required to be in the up position (open), even if the vacuum manifold is at atmospheric pressure. This is to be consistent with maintenance procedures that rely on the inlet valve lever being closed at the appropriate step to prevent the accidental loss of vacuum. If the vacuum is accidentally lost the system may be damaged. At a minimum, the components that were under vacuum might have to be cleaned. ▲



Labeled components: 1=inlet valve plug, 2=vacuum manifold, 3=inlet valve lever, 4=guide bar, 5=entry housing, 6=inlet valve knob, 7=inlet valve block, 8=foreline hose connection, 9=12 pin feedthrough, 10=vacuum manifold probe plate, 11=ball valve housing, 12= inlet valve solenoid

Figure 3-27. Inlet valve components (ion volume tool not shown)

- b. Remove all connectors between the components on the vacuum manifold probe plate (item #10, [Figure 3-27](#) on [page 3-34](#)) and the ETD Control PCB ([Figure 1-16](#) on [page 1-22](#)).
- c. Remove the valve shield from the vacuum manifold probe plate ([Figure 3-28](#)) by loosening the four screws at the corners of the shield.

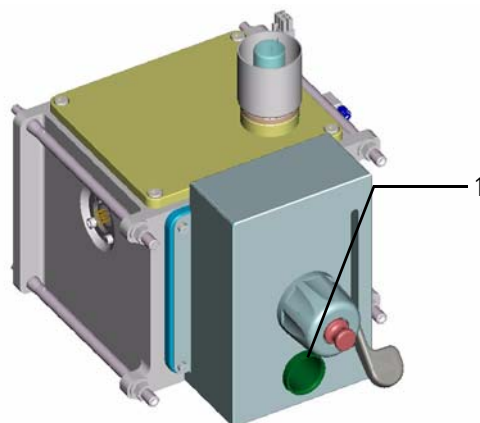


Figure 3-28. Valve shield (1) covering the vacuum manifold probe plate

- d. Remove the foreline hose on the source from its connection (Figure 3-29 and item 8 in Figure 3-27).



Figure 3-29. Removing the foreline hose from its connection

- e. Remove the four screws holding the vacuum manifold probe plate (Figure 3-30 and item 10 in Figure 3-27 on page 3-34). Support the plate with your hand as shown in Figure 3-30. Arrows point to the four hex screw locations (items 1–4).

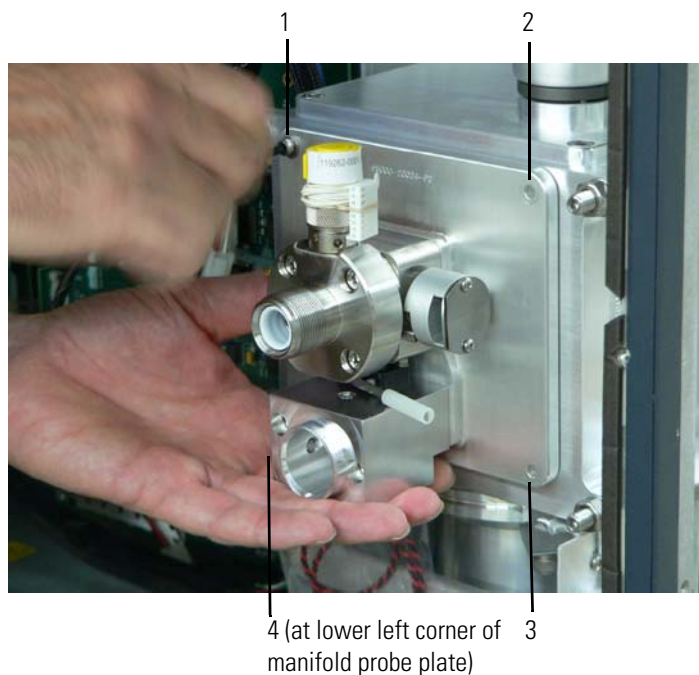


Figure 3-30. Unscrewing the vacuum manifold probe plate

- f. Remove the vacuum manifold probe plate ([Figure 3-31](#) on [page 3-36](#)).

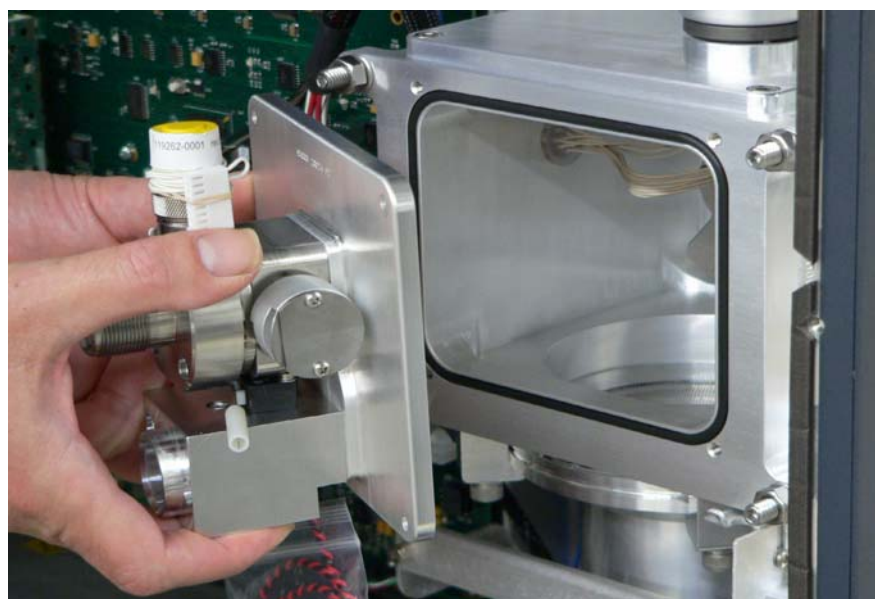
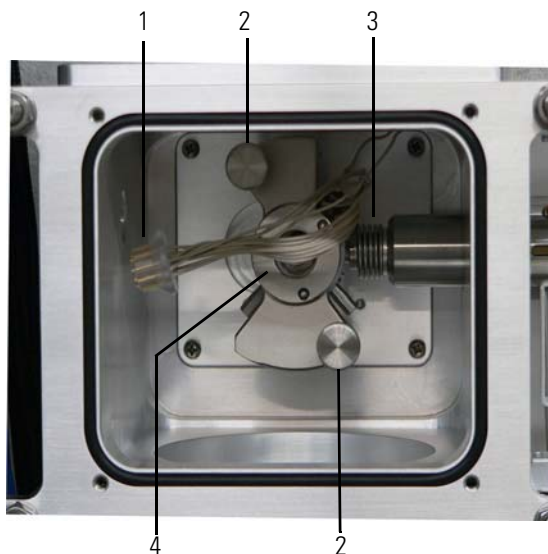


Figure 3-31. Removing the vacuum manifold probe plate

- g. Unplug the 12-pin feedthrough harness from the feedthrough ([Figure 3-32](#)).



Labeled components: 1=unplugged 12 pin feedthrough, 2=thumbscrews, 3=transfer line bellows, 4=ion source assembly

Figure 3-32. Interior of vacuum manifold

- h. Remove the ion source assembly from the vacuum manifold (Figure 3-33 on page 3-38) by first loosening the ion source thumbscrews (item 2 in Figure 3-32).
- i. Second, as you remove the ion source assembly (item 1 in Figure 3-33) gently shift it to the left (arrow 2 in Figure 3-33) before and while pulling it out. This will allow the ion source assembly to disengage from the transfer line bellows (item 3 in Figure 3-33) as it is removed. Alternatively, gently depress the transfer line bellows (Figure 3-32 on page 3-37) to disengage it from the ion source assembly.

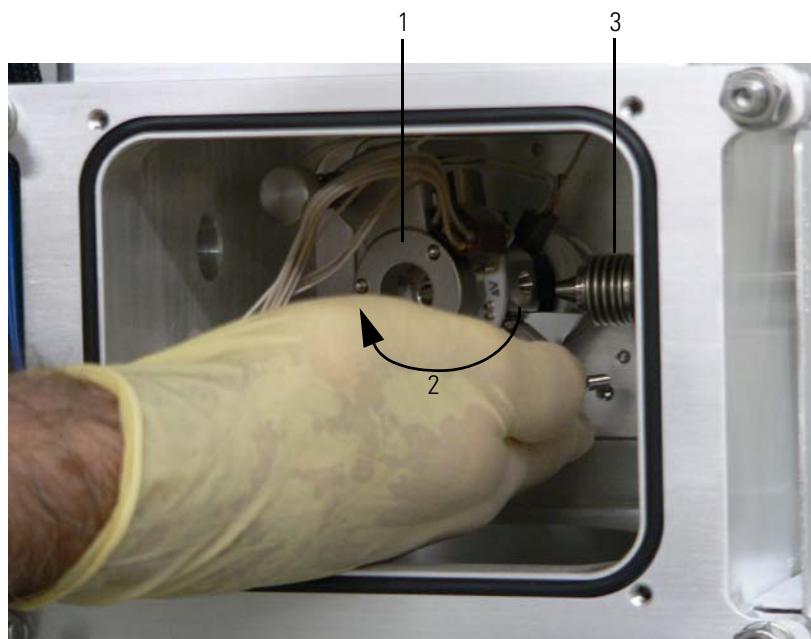
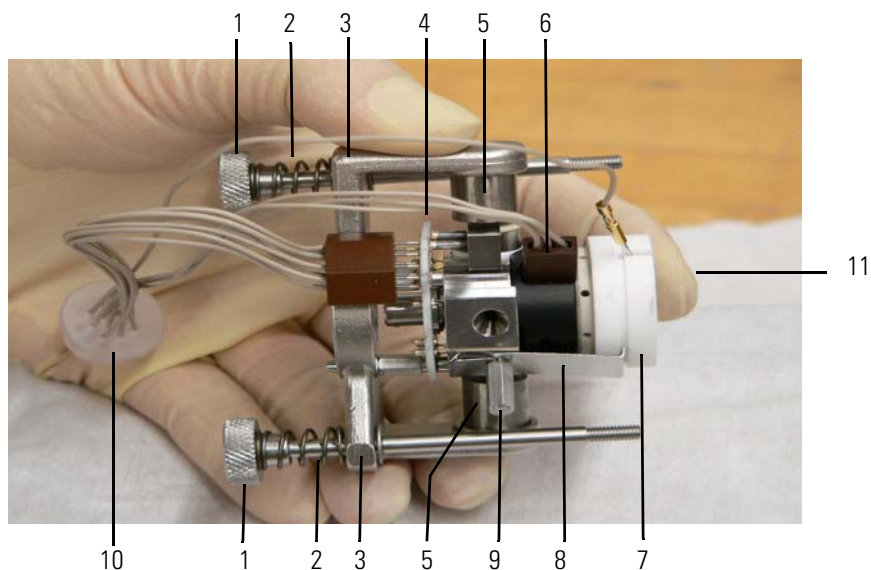


Figure 3-33. Removing the ion source assembly from the vacuum manifold^a

^aThe ion source assembly (item 1) is gently shifted to the left (arrow 2) to allow the ion source assembly to disengage from the transfer line bellows (item 3) as it is removed.

The ion source assembly is held together with a clip (item 8 in [Figure 3-34](#) on [page 3-39](#)). However, it is necessary to keep the tips of your gloved fingers on both the front edge of the ceramic lens holder (item 11 in [Figure 3-34](#)) and the back of the magnet yoke (item 3 in [Figure 3-34](#)) when you handle the ion source assembly. This prevents unsecured components inside of the ceramic lens holder from falling out.

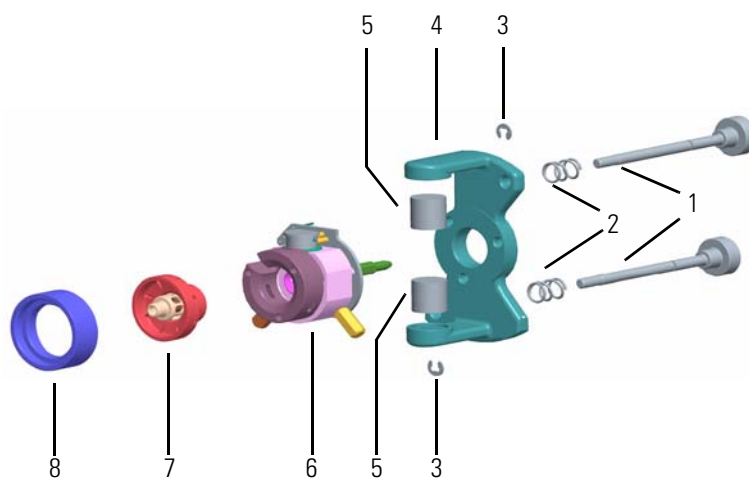
Caution When handling the ion source assembly, it is important to handle it with gentle finger pressure on each end (as instructed above) to keep unsecured components from falling out of the assembly. ▲



Labeled components: 1=thumbscrew, 2=springs, 3=magnet yoke, 4=Ion Source PCB, 5=magnets, 6=ion source block, 7=ceramic lens holder, 8=spring clip, 9=spring clip thumb screw, 10=12 pin feedthrough harness, 11=finger over front edge of ceramic lens holder to keep unsecured components from falling out of the ion source assembly.

Figure 3-34. Ion source assembly

An exploded view of the ion source assembly is shown in [Figure 3-35](#).



Labeled components: 1=thumbscrews, 2=springs, 3=E-clips, 4=magnet yoke, 5=magnets, 6=ion source, 7=ion source lens assembly, 8=ceramic lens holder

Figure 3-35. Ion source assembly exploded view

3. Separate the magnet yoke and the ion source.
4. Remove the ion source lens assembly from the ceramic lens holder ([Figure 3-35](#)).

5. Clean the ion source lens assembly according to the procedure in “[Cleaning Stainless Steel Parts](#)” on [page 3-14](#). Pay particular attention to the areas inside the tube and around the holes in the lens assembly.
6. Replace the ion source assembly:
 - a. Insert the lens assembly into the ceramic lens holder.
 - b. Reassemble the ion source assembly.
 - c. Reinstall the ion source assembly into the vacuum manifold by following [step 2](#) in reverse order.
7. Restore the ETD Module to operational status. See “[Starting Up the System after a Shutdown](#)” on [page 2-9](#).

Cleaning the Ion Source Block

If cleaning the ion volume and ion source lens assembly does not restore system performance, you might need to clean the ion source block. Generally, you need to clean the ion source block no more than once every six months.

Supplies needed for cleaning the ion source:

- Cleaning supplies
- Gloves (clean, lint-free, and powder-free)
- Lint-free cloth

❖ To clean the ion source block

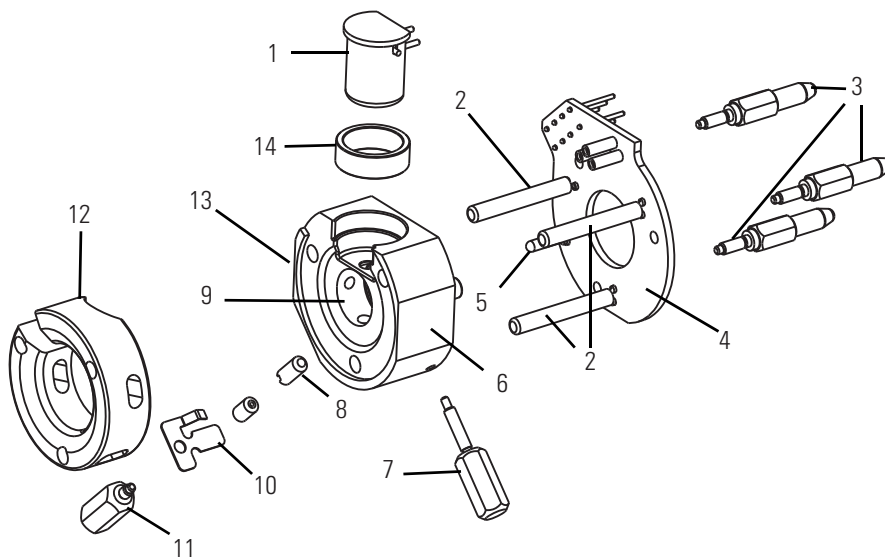
1. Prepare the ETD Module for maintenance:
 - a. Prepare a clean work area by covering the area with lint-free cloth.
 - b. Shut down and vent the Orbitrap Velos Pro ETD mass spectrometer. (See “[Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely](#)” on [page 2-7](#).)

Caution Shut down and unplug the Orbitrap Velos Pro ETD mass spectrometer before proceeding with the next steps of this procedure. ▲

- c. Remove the ion source assembly by following the procedures in [step 2](#) in “[Cleaning the Ion Source Lens Assembly](#)” on [page 3-33](#).

Note Wear clean, lint- and powder-free gloves when you handle parts inside the vacuum manifold. ▲

2. Disassemble the ion source assembly (Figure 3-34 on page 3-39 and Figure 3-35 on page 3-39), remove and disassemble the ion source (Figure 3-36).
 - a. Remove the magnet yoke and the ion source block with the ion source lens assembly.
 - b. Remove the ion source lens assembly.
 - c. Remove the ion source.



Labeled components: 1=ion source filament, 2=cartridge heaters, 3=base studs (3×), 4=Ion Source PCB, 5=temperature sensor, 6=ion source block, 7=ion volume key thumbscrew, 8=ion volume pin, 9=ion volume, 10=spring clip, 11=spring clip thumb screw, 12=heater ring, 13=sample inlet aperture (in side of item 6), 14=ceramic spacer

Figure 3-36. Ion source, exploded view

- d. Remove the three base studs (item 3 in Figure 3-36). Be careful not to damage the leads on the Ion Source PCB (item 4 in Figure 3-36).
- e. Gently remove the Ion Source PCB (item 4 in Figure 3-36) from the ion source by loosening the spring clip thumbscrew (item 11 in Figure 3-36) and the spring clip (item 10 in Figure 3-36) and sliding the three cartridge heaters and the temperature sensor (items 2 and 5 in Figure 3-36) off the ion source and pulling the filament (item 1 in Figure 3-36) straight away from the three filament connectors on the Ion Source PCB. Do not bend or twist the cartridge heaters or temperature sensor.
- f. Remove the filament and ceramic spacer (items 15 and 1 in Figure 3-36) from the ion source block (item 6 in Figure 3-36).

- g. Remove the ion volume key thumbscrew (item 7 in [Figure 3-36](#) on [page 3-41](#)).

Note It is not necessary to remove the ion volume pin (item 8 in [Figure 3-36](#)). If you remove it, you should reinsert it just far enough so the ball will keep an ion volume (item 9 in [Figure 3-36](#)) from falling out. If the ball extends too far, the ion volume will be difficult to remove. ▲

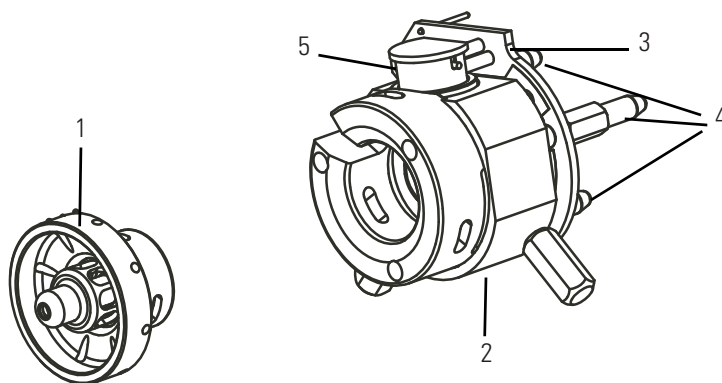
3. Clean the ion source parts and replace the ion source assembly:
 - a. Clean each component of the ion source, as described in “[Cleaning Stainless Steel Parts](#)” on [page 3-14](#) and “[Cleaning Non-Stainless Steel or Hybrid Parts](#)” on [page 3-17](#).
 - b. Reassemble the ion source block.
 - c. Reassemble the ion source assembly.
 - d. Reinstall the ion source assembly into the vacuum manifold by following [step 2](#) of “[Cleaning the Ion Source Lens Assembly](#)” on [page 3-33](#) in reverse order.
4. Restore the ETD Module to operational status. See “[Starting Up the System after a Shutdown](#)” on [page 2-9](#).

Replacing the Ion Source Filament

The number of ions produced in the ion source is approximately proportional to the filament emission current. If you notice that ion production is low, this might indicate that the filament has failed and needs to be replaced. If the measured emission current is substantially less than the value that the emission current is set to, or if the measured emission current is decreasing over time, then the filament has failed or is failing and needs to be replaced. The ion source filament assembly is shown in [Figure 3-37](#) on [page 3-43](#).

Supplies needed for replacing the ion source filament:

- Filament Assembly DSQ II (P/N 120320-0030)
- Gloves, clean, lint-free, and powder-free
- Protective eyewear
- Lint-free cloth
- Forceps or dental pick



Labeled components: 1=ion source lens assembly, 2=ion source block, 3=Ion Source PCB, 4=base studs (3x), 5=ion source filament

Figure 3-37. Ion source lens assembly and ion source

❖ **To replace the ion source filament**

1. Prepare the ETD Module for maintenance.
 - a. Prepare a clean work area by covering the area with lint-free cloth.
 - b. Shut down and vent the Orbitrap Velos Pro ETD mass spectrometer. (See “[Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely](#)” on page 2-7.)

Caution Shut down and unplug the Orbitrap Velos Pro ETD mass spectrometer before proceeding with the next steps of this procedure. ▲

- c. Remove the ion source assembly by following the procedures in step 2 in “[Cleaning the Ion Source Lens Assembly](#)” on page 3-33.

Note Wear clean, lint- and powder-free gloves when you handle parts inside the vacuum manifold. ▲

2. Disassemble the ion source assembly ([Figure 3-34](#) on page 3-39 and [Figure 3-35](#) on page 3-39), remove and disassemble the ion source ([Figure 3-36](#) on page 3-41).
 - a. Remove the ion source lens assembly (item 1 in [Figure 3-37](#)).
 - b. Remove the three base-studs (item 3 in [Figure 3-36](#), item 4 in [Figure 3-37](#)).
 - c. Remove the filament assembly (items 1 and 15 in [Figure 3-36](#), item 5 in [Figure 3-37](#)) and ion source block (item 2 in [Figure 3-37](#)) from the three filament connectors and cartridge heaters (item 2 in [Figure 3-36](#)) on the Ion Source PCB (item 4

in [Figure 3-36](#)) according to the procedure in [step e](#) of “Cleaning the Ion Source Block” on [page 3-41](#) .

Note Now is a good time to clean the ion volume and ion source lenses. ▲

3. Inspect and install a new filament assembly:
 - a. Turn the filament assembly over and, using a strong light and a magnifying glass, verify that the filament wire is centered in the electron lens hole. If necessary, carefully use forceps (or a dental pick) to adjust the filament wire. [Figure 3-38](#) shows the centered filament wire as seen from the bottom of the filament through the electron lens hole.

Note A bent filament can lead to a low or absent anion signal. If the filament wire is bent, not centered, or otherwise damaged, you must replace the filament assembly. ▲

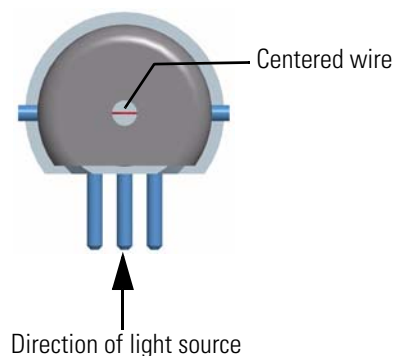


Figure 3-38. Filament wire as seen from the bottom of the filament through the electron lens hole

- b. Insert the filament into the ceramic spacer of the ion source block (item 14 in [Figure 3-36](#) on [page 3-41](#)).
 - c. Align the filament leads with the Ion Source PCB connectors and gently press the leads into the connectors. Normally, there is a small gap (about 0.020 in or 0.50 mm) between the filament and the connectors. The gap allows the ceramic filament centering ring (spacer) to properly position and align the electron lens hole with the ion volume.
 - d. Reinstall the three base-studs (item 3 in [Figure 3-36](#), item 4 in [Figure 3-37](#) on [page 3-43](#)).
4. Reassemble ion source and ion source assembly.
5. Insert the ion source assembly into the vacuum manifold.

6. Restore the ETD Module to operational status. See [“Starting Up the System after a Shutdown”](#) on [page 2-9](#).

Note Tune Plus provides an evaluation procedure for CI gas pressure under **Diagnostics > Diagnostics > Tools > System evaluation > Reagent CI gas pressure evaluation**. Thermo Fisher Scientific recommends performing this procedure after replacing the filament and/or the ion volume. ▲

Replacing Inlet Valve Components

This topic provides the procedure for replacing inlet valve components. Perform this procedure when the inlet valve seal or the inlet valve is being replaced.

Tools and supplies needed for replacing inlet valve components:

- Inlet Valve Seal Kit (P/N 119265-0003)
- Lint-free cloth
- Wrench, open-ended, 5/16-in
- Wrench, Allen, 4 mm

❖ To replace inlet valve components

1. Pull down the inlet valve lever to close the inlet valve. See [Figure 3-39](#).

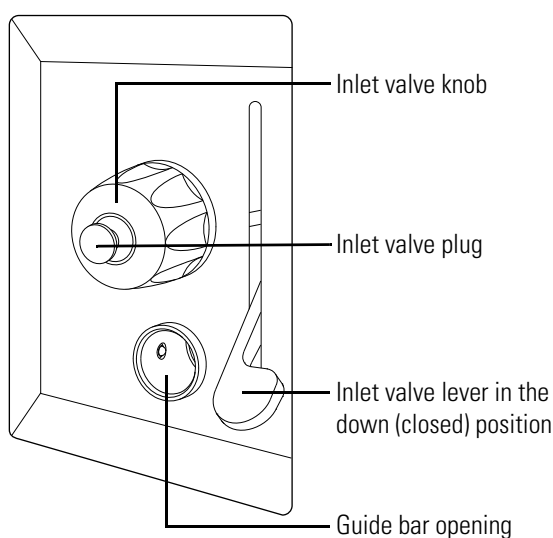


Figure 3-39. Inlet valve components

2. Rotate the inlet valve knob counter-clockwise until you can easily remove the inlet valve plug.

3. Continue to rotate the valve knob until you can remove it.

The inlet valve knob has a stainless-steel ferrule inside. Keep the inlet valve knob and ferrule together.

4. Pull out the knob on the inlet valve seal tool (P/N 119283-0001)¹, so that the knob is loose. See [Figure 3-40](#).

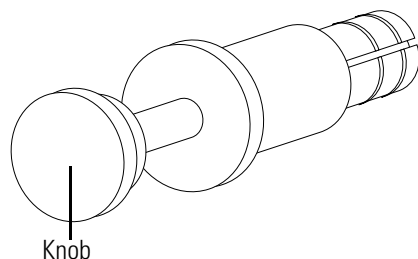


Figure 3-40. Inlet valve seal tool

5. Insert the tool straight into the inlet valve. See [Figure 3-41](#).

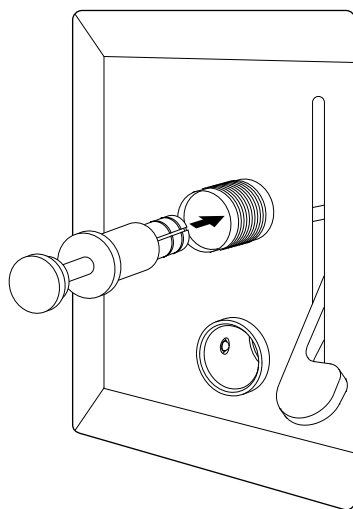


Figure 3-41. Inlet valve seal tool inserted in the inlet valve

Caution Do not scratch the surface of the seal. Use only the inlet valve seal tool to remove or install an inlet valve seal. ▲

6. Press in the knob on the tool until it stops.
7. Remove the tool. The inlet valve seal should be on the tool. See [Figure 3-42](#) on [page 3-47](#).

¹Item contained in Inlet Valve Seal Kit (P/N 119265-0003).

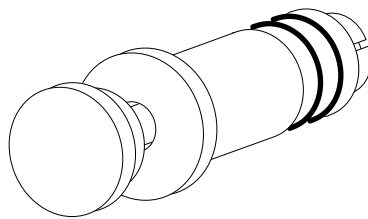


Figure 3-42. Inlet valve seal on the inlet valve seal tool

8. Loosen the knob to disengage the seal. See [Figure 3-43](#).

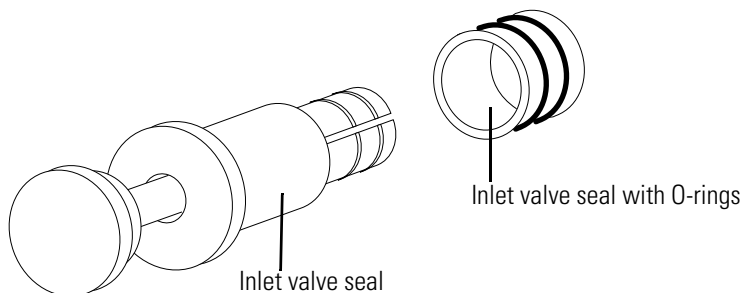


Figure 3-43. Inlet valve seal disengaged from tool

9. Discard the seal with its O-rings in place.
10. Place one O-ring (P/N 3814-6530)¹ into each of the two slots on the inlet valve seal (P/N 119683-0100)¹.
11. Place the new inlet valve seal onto the inlet valve seal tool.
12. Insert the inlet valve seal tool into the inlet valve into it stops.
13. Remove the tool. The O-rings on the valve seal secure the inlet valve seal in the opening.
14. Reinstall the ferrule, knob, and plug into the inlet valve opening.

¹Item contained in Inlet Valve Seal Kit (P/N 119265-0003).

Replacing the Reagent Vials

A significant drop of the m/z 202 signal within one hour with the mission current at the correct level indicates a low reagent supply. In this case, you should replace the fluoranthene vial.

Changing the reagent vials requires that the Orbitrap Velos Pro ETD mass spectrometer be placed in Service mode after the vials have cooled. (Vial cooling is done in Off condition.) See the following sections for procedures to be used to change the reagent vials:

- “[Placing the Instrument in Off Condition and Service Mode](#)” on page 3-48
- “[Installing/Exchanging the Reagent Vials](#)” on page 3-51

The ETD reagent vials are designed to keep the ETD reagent out the lab environment. Removal and replacement of the ETD reagent vials when they are not empty causes excessive puncturing of the septums and reduces their integrity. This could result in ETD reagent (fluoranthene) entering the lab environment. Prevent this from occurring by removing and replacing the ETD reagent vials only when they are empty.

Caution To preserve the integrity of the ETD reagent vial septums, remove and replace the ETD reagent vials only when they are empty. Do not reinstall used vials. ▲

Note Store and handle all chemicals in accordance with standard safety procedures. The Material Safety Data Sheet (MSDS) describing the chemicals being used should be freely available to lab personnel for them to examine at any time. Material Safety Data Sheets (MSDSs) provide summarized information on the hazard and toxicity of specific chemical compounds.

MSDSs also provide information on the proper handling of compounds, first aid for accidental exposure, and procedures for cleaning spills or dealing with leaks. Producers and suppliers of chemical compounds are required by law to provide their customers with the most current health and safety information in the form of an MSDS. Read the MSDS for each chemical you use. Dispose of all laboratory reagents in the appropriate manner (see the MSDS). ▲

Safety information about fluoranthene is given in [Appendix A: “Fluoranthene”](#).

Placing the Instrument in Off Condition and Service Mode

The power switches control power to the Orbitrap Velos Pro ETD mass spectrometer (MS and ETD Module). The ETD Module power switches control the power to the ETD Module only. When the

Orbitrap Velos Pro ETD MS is fully operational (all systems On), the MS Main Power switch is in the On position and the FT Electronics switch is in the operating (On) position.

Normally, the ETD Module Power and Service switches remain On. Use the FT Electronics switch on the MS unit to place the Orbitrap Velos Pro ETD mass spectrometer in Service mode. Turn On and Off the instrument (both ETD Module and MS) with the MS Main Power switch.



Warning Burn Hazard. When mass spectrometer and ETD Module system are turned On, the flow restrictor, the transfer line heaters, and the ion source heater can be at 160 °C. The vial heaters can be at 108 °C (or set point). Do not attempt to replace reagent vials or to service heated components until you have determined that they have cooled to a safe temperature for handling. ▲

Note The instructions that follow assume that no analyte is flowing into the API source. ▲

❖ **To place the Orbitrap Velos Pro ETD mass spectrometer in Off Condition and Service mode and to verify that the vials are safe to handle**

1. If the Tune Plus window is not already open, choose **Start > Programs > Thermo Instruments > LTQ > LTQ Tune** from the taskbar. The Tune Plus window appears. (See [Figure 3-10](#) on [page 3-22](#).)



On

Off

Standby

You can determine the state of the MS detector by observing the state of the On/Standby button on the Control/Scan Mode toolbar. The three different states of the On/Standby button are shown at the left.

2. Choose **Control > Off** from the Tune Plus pull-down menu to place the system in Off condition. When the mass spectrometer is in Off condition, the Orbitrap Velos Pro ETD MS turns off the ion source sheath gas, auxiliary gas, high voltage, and all of the ETD Module heaters.

Note It is important to choose **Control > Off** from the Tune Plus pull-down menu in order to shut down all of the ETD Module heaters. ▲



3. Click the reagent ion source portion of the instrument control icon at the top of the Tune Plus window. The Reagent Ion Source dialog box appears ([Figure 3-44](#)).

Observe the temperature of Vial 1 in the Actual column of the Reagent Ion Source dialog box (Figure 3-44). Nitrogen cooling gas will flow until the vial reaches 70 °C. (See “Turning Off the Reagent Ion Source: What to Expect” on page 2-14.) Allow up to 90 minutes for the vial temperature to reach ambient temperature (about 30 °C).

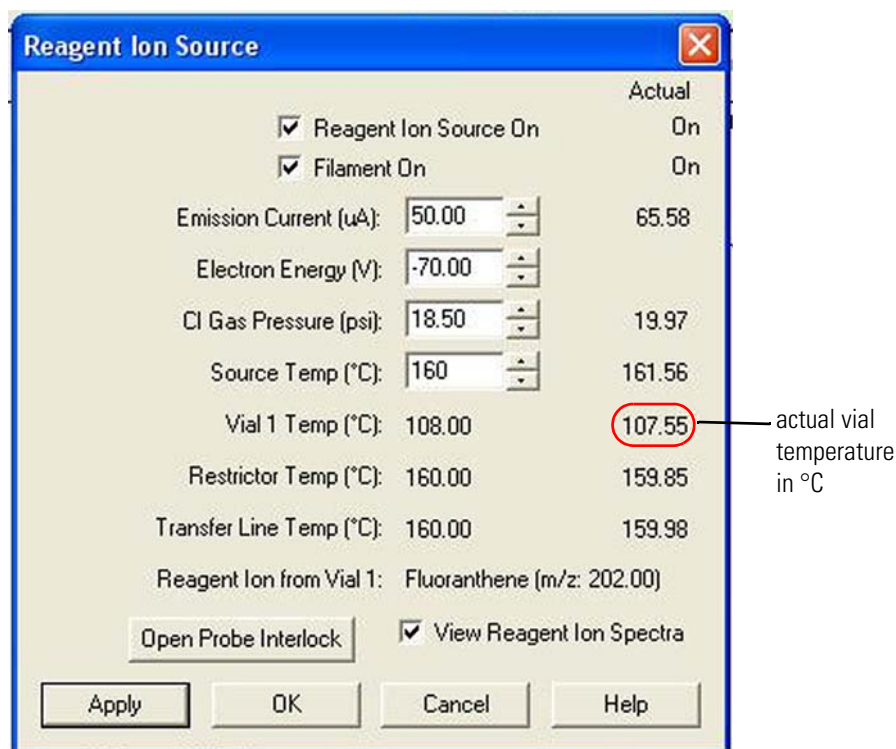


Figure 3-44. Reagent Ion Source dialog box



Warning Burn Hazard. Do not attempt to handle the vials or vial holders when the cooling nitrogen stops. They are still too hot to handle when the cooling nitrogen stops at a vial temperature of 70 °C. Allow the vials to cool to about 30 °C (allow up to 90 minutes after the cooling gas stops) before proceeding with the next step and handling the vials. ▲

4. Toggle the FT Electronics switch to Service mode (Off) when the vial has reached a temperature that is safe for handling (about 30 °C). Toggling the FT Electronics switch to Service mode turns off all components except the turbomolecular pumps and the forepumps in both the mass spectrometer and the ETD Module.

Note Do not place the ETD Module Service switch into its Service mode (Off) position while the MS switches are left in their On positions. This could cause communication problems between the MS and the ETD Module. The ability to control the Service mode for both the MS and the ETD Module at one point (at the FT Electronics switch) is a safety feature. ▲



Warning Burn Hazard. Do not place the system in Service mode until the vials reach a safe temperature (about 30 °C). System temperature monitoring will stop when the system is placed in Service mode. Do not attempt to handle the vials, the vial holders, or the heater assembly until a safe temperature is reached (about 30 °C). ▲

The Orbitrap Velos Pro ETD mass spectrometer is now in Service mode and the vials are at a safe temperature for handling.

Installing/Exchanging the Reagent Vials

After the reagent vial heaters have cooled to room temperature, the reagent vials are ready to be installed or exchanged.

❖ To install or exchange the reagent vials

1. Remove the back panel from the ETD Module. (See “[Removing the ETD Main Access Panel](#)” on [page 3-18](#).) This exposes the reagent inlet source heating unit, which has its own cover ([Figure 3-45](#) on [page 3-52](#)).



Warning Burn Hazard. Follow the procedures described in “[Placing the Instrument in Off Condition and Service Mode](#)” on [page 3-48](#) before removing the back panel of the ETD Module. Removing the back panel before the system is placed in Service mode will open the panel electrical interlocks and stop all system activity including temperature monitoring. In the absence of temperature monitoring, you might attempt to handle the vials before it is safe to do so. ▲

2. Make sure that the vial heater cover is cool to the touch.



Warning Burn Hazard. The vial heaters can be at 108 °C (or set point). Allow sufficient time for the vials to cool (up to 90 minutes) and then place the system in Service mode. (See “[Placing the Instrument in Off Condition and Service Mode](#)” on [page 3-48](#).) Verify that the vial heater cover is safe to handle before attempting to remove the vial holders and reagent vials. ▲

User Maintenance

Maintenance of the ETD Module

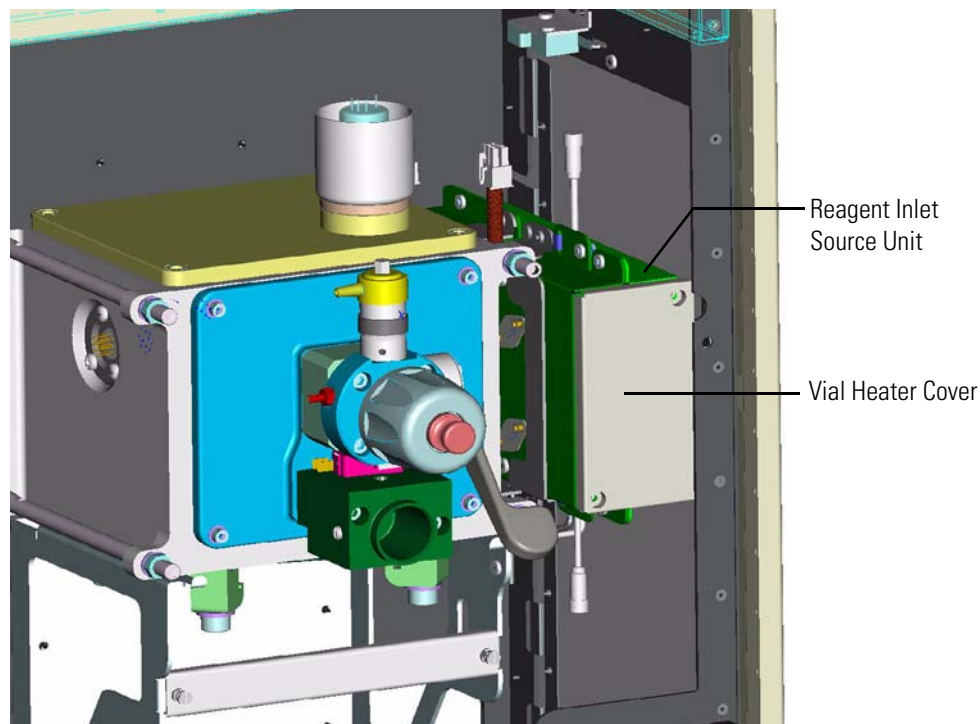
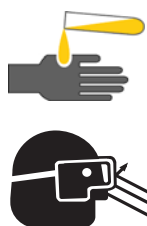


Figure 3-45. ETD Module with back panel removed

3. Put on a pair of new, white nitrile clean room gloves and protective eye wear.
4. Using a Phillips screwdriver, remove the screws in the vial heater cover. The vial heater cover is located on the right side of the ETD Module as you view it from the back of the Orbitrap Velos Pro ETD mass spectrometer (Figure 3-45).
5. Remove the vial holder by gently pulling it out of the vial heater.
6. Remove the empty vial if it is present. The vial holder is a cylindrical tube with a handling knob at one end and ribs along its length. See Figure 3-46. These ribs prevent the vial holder from rotating once it is placed into the vial heater. Figure 3-47 on page 3-53 shows the tab and ribs of a vial holder in the vial heater.

Note Dispose of an empty fluoranthene vial in accordance with its MSDS. ▲



Warning Avoid exposure to potentially harmful materials.

Always wear protective gloves and safety glasses when you handle solvents or corrosives. Also contain waste streams and use proper ventilation. Refer to your supplier's Material Safety Data Sheet (MSDS) for proper handling of a particular compound. ▲



Figure 3-46. Reagent vials with holders

7. Take a vial containing the ETD reagent (fluoranthene) from its box and remove the aluminum tab from the top of the vial's crimp seal.
8. Put the vial into a vial holder.

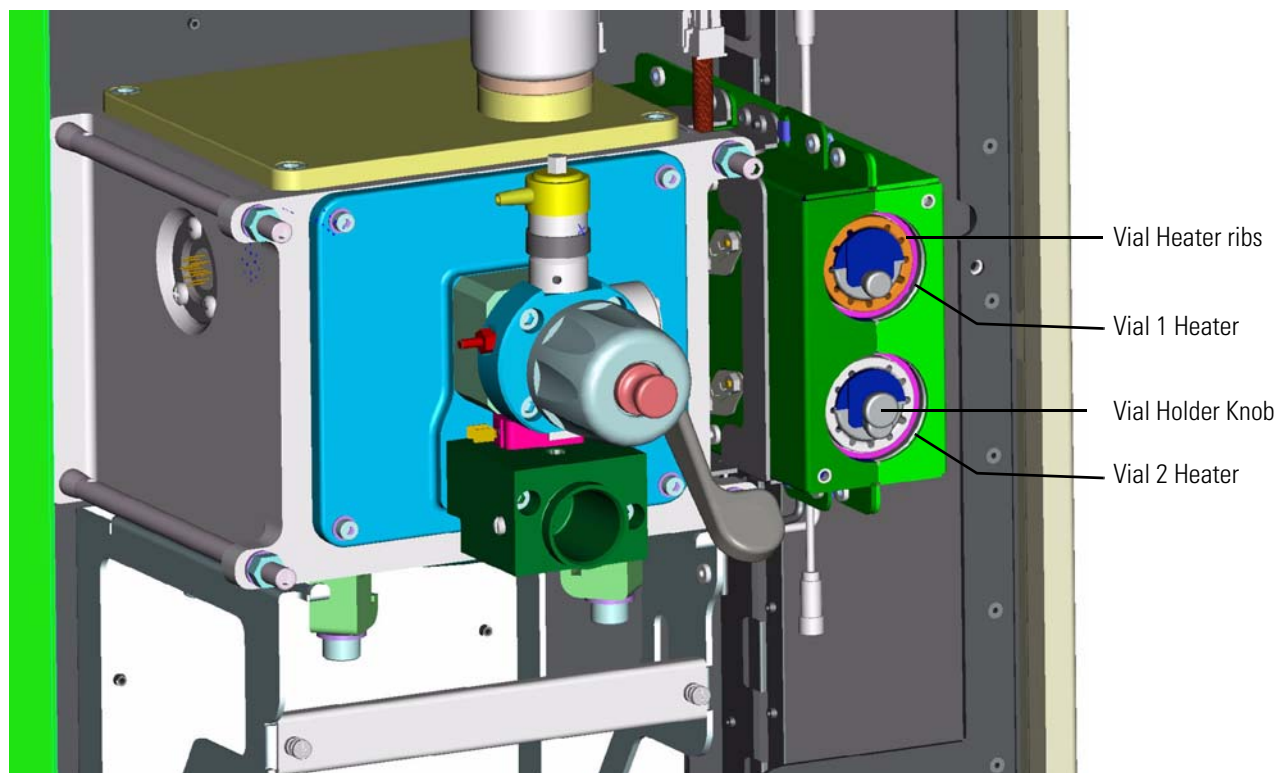


Figure 3-47. ETD Module with vial heater cover removed

9. Place this ETD reagent vial and its vial holder into the Vial 1 heater (top vial heater). Gently slide the vial holder into the vial heater.
10. Place the empty vial from the box into the other vial holder if an empty vial is not already installed.
11. Place this empty vial and its vial holder into the Vial 2 heater (bottom vial heater) if an empty vial is not already installed.



Warning Health Hazard. The empty vial in the Vial 2 heater is an integral part of the carrier/CI gas system. It is necessary to keep the carrier/CI gas system closed to the laboratory. If no vial is placed in the Vial 2 heater, the carrier/CI gas containing the reagent might escape to the laboratory causing a safety problem. ▲

Caution If no vial is placed in the Vial 2 heater, the ETD Module will not operate correctly and the filament will burn out. ▲

12. Reinstall the vial heater cover over the vial heaters.
13. Reinstall the back panel of the ETD Module. See “[Removing the ETD Main Access Panel](#)” on [page 3-18](#). The ETD Module will not turn on unless the back panel is installed.
14. Start the system:
 - a. Toggle the FT Electronics switch to the On position. The system will boot to Standby mode. Then ion source heater, flow restrictor, and transfer line heaters will start heating. Monitor these temperatures in the Status View on the right side of the Tune Plus window. (See [Figure 3-10](#) on [page 3-22](#).) They will have green check marks when they have reached their operating temperatures.
 - b. Select the Reagent Ion Source On check box in the Reagent Ion Source dialog box when the ion source heater, flow restrictor, and transfer line heaters are at their operating temperatures. (See [Figure 3-44](#) on [page 3-50](#).)

The Orbitrap Velos Pro ETD mass spectrometer is now ready for use.

Changing the Reagent Ion Source Flow Restrictors

❖ To change the reagent ion source flow restrictors

1. Shut down completely the instrument according to the procedures in “[Shutting Down the Orbitrap Velos Pro Mass Spectrometer Completely](#)” on [page 2-7](#).



Warning Burn Hazard. The reagent vial heaters can be 108 °C (or set point), the flow restrictor, the transfer line heaters, and the ion source heater can be at 160 °C. These components may be too hot to touch. Verify that all of these components are safe to touch before handling them. ▲

2. Remove the ETD side access panel according to the instructions in “Removing the ETD Side Access Panel” on page 3-19.

Caution The ETD side access panel is interlocked with the ETD Module power. When the ETD side access panel is removed, all power to the ETD Module will be turned off. ▲

3. Remove the four screws that hold the reagent inlet cover in place.



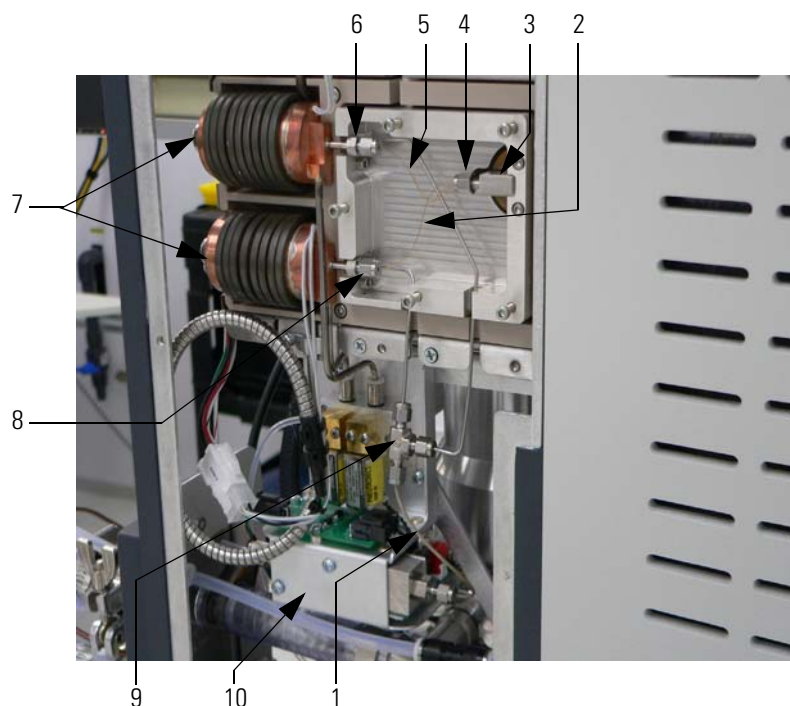
Warning Burn Hazard. The reagent inlet cover may be too hot to touch. Verify that the reagent inlet cover is at or near room temperature before handling it. ▲

4. Remove the five screws that hold the restrictor oven cover in place.



Warning Burn Hazard. The restrictor oven cover may be too hot to touch. Verify that the restrictor oven cover is at or near room temperature before handling it. ▲

5. Replace both fused silica restrictors and their ferrules by removing their Swagelok® nuts. See [Figure 3-48](#).



Labeled components: 1=PEEKsil[®] tubing, 2=fused silica tubing from lower oven, 3=transfer line inlet, 4=Swagelok fitting with two hole ferrule, 5=fused silica tubing from upper oven, 6=upper oven Swagelok fitting with a single hole ferrule, 7=vial 1 (upper) and vial 2 (lower) heaters, 8=lower oven Swagelok fitting with a single hole ferrule, 9=Tee below reagent inlet assembly, 10=gas valves

Figure 3-48. Reagent inlet assembly

6. Thread two pieces of fused silica tubing (P/N 98000-20060) into the two hole ferrule (P/N 00101-08-00006) from the Installation Kit for the Reagent Inlet Module (P/N 98000-62006). Place a Swagelok fitting over the ferrule.
7. Insert the two hole ferrule and Swagelok fitting from [step 6](#) onto the transfer line inlet (item 3 in [Figure 3-48](#)) and tighten the Swagelok fitting.
8. Thread the opposite end of one of the pieces of fused silica tubing into a single hole ferrule. Place a Swagelok fitting over the ferrule.
9. Insert the ferrule and Swagelok fitting from [step 8](#) on to one of the oven outlets (items 6 or 8 in [Figure 3-48](#)) and tighten the Swagelok fitting.

Caution Do not overtighten the ferrules. The ferrules may loosen after they are first heated. If this occurs retighten them if necessary. ▲

10. Repeat [step 9](#) for the other oven.

11. Replace the restrictor oven cover and reagent inlet cover removed in [step 3](#) and [step 4](#).
12. Loosen the Swagelok fitting connecting the PEEKsil® tubing (item 1 in [Figure 3-48](#)) to the Tee below the reagent inlet assembly (item 9 in [Figure 3-48](#)) and from the gas valves (item 10 in [Figure 3-48](#)).
13. Replace the old PEEKsil tubing with new PEEKsil tubing (P/N 00109-02-00020) from the Installation Kit for the Reagent Inlet Module (P/N 98000-62006).
14. Close the ETD Module and restart the instrument. Follow the procedures given in [“Starting Up the System after a Shutdown”](#) on [page 2-9](#).

Cleaning the Fan Filters of the ETD Module

You need to clean the fan filters every four months. The fan filters are located at the rear of the ETD Module on the left side (as viewed from the back of the ETD Module). See [Figure 3-49](#).

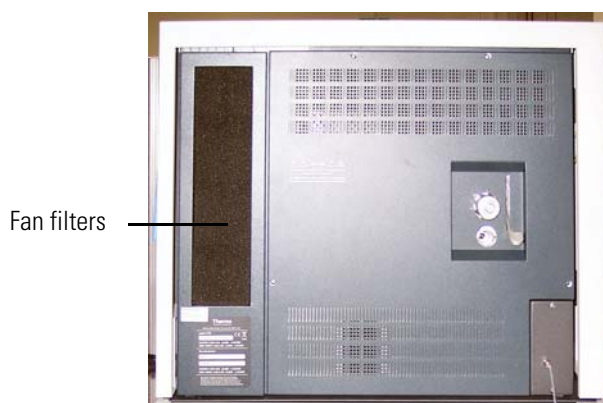


Figure 3-49. ETD Module, top panel

❖ To clean the fan filters of the ETD Module

1. Remove the fan filter from the rear of the ETD Module by pulling it out of the fan filter bracket.
2. Wash the fan filters in a solution of soap and water.
3. Rinse the fan filters with tap water.
4. Squeeze the water from the fan filters and allow them to air dry.
5. Reinstall the fan filter in the fan filter bracket.

Maintenance of the Cooling Circuit

The recirculating chiller and the water filter require maintenance on a regular basis.

Maintenance for the Recirculating Chiller

For the NESLAB ThermoFlex™ 900 recirculating chiller, the checks described in this section should be carried out on a regular basis.

Note For further information and maintenance instructions, refer to the manufacturer's manual supplied with the instrument. ▲

Reservoir

Periodically inspect the fluid inside the reservoir. If cleaning is necessary, flush the reservoir with a cleaning fluid compatible with the circulating system and the cooling fluid.

The cooling fluid should be replaced periodically. Replacement frequency depends on the operating environment and amount of usage.



Warning Burn Hazard. Before changing the operating fluid, make sure it is at safe handling temperature. ▲

Fluid Bag Filter

The ThermoFlex 900 recirculating chiller installed in the cooling circuit of the instrument is equipped with a fluid bag filter, which needs to be replaced on a regular basis. Replacement bags are available from Thermo Fisher Scientific.

Condenser Filter

To prevent a loss of cooling capacity and a premature failure of the cooling system, clean the condenser filter regularly. If necessary, replace it.

Replacing the Water Filter Cartridge

The filter removes particulate matter in the cooling system that might damage the flow sensor. The filter cartridge should be replaced annually or as necessary. A replacement is available from Thermo Fisher Scientific (P/N 1284050).

❖ **To replace the water filter cartridge**

1. Place the system in Standby condition as described on [page 2-4](#).
2. Turn off the chiller.
3. The filter is installed on the left instrument side, in the cooling lines of the Orbitrap system between the Peltier element and the flow sensor. See [Figure 3-50](#). Quick couplings connect the filter assembly to the hoses of the cooling lines. Press the thumb latch of each quick coupling to release it; valves in the couplings prevent the water from leaking out. Then remove the complete filter assembly.

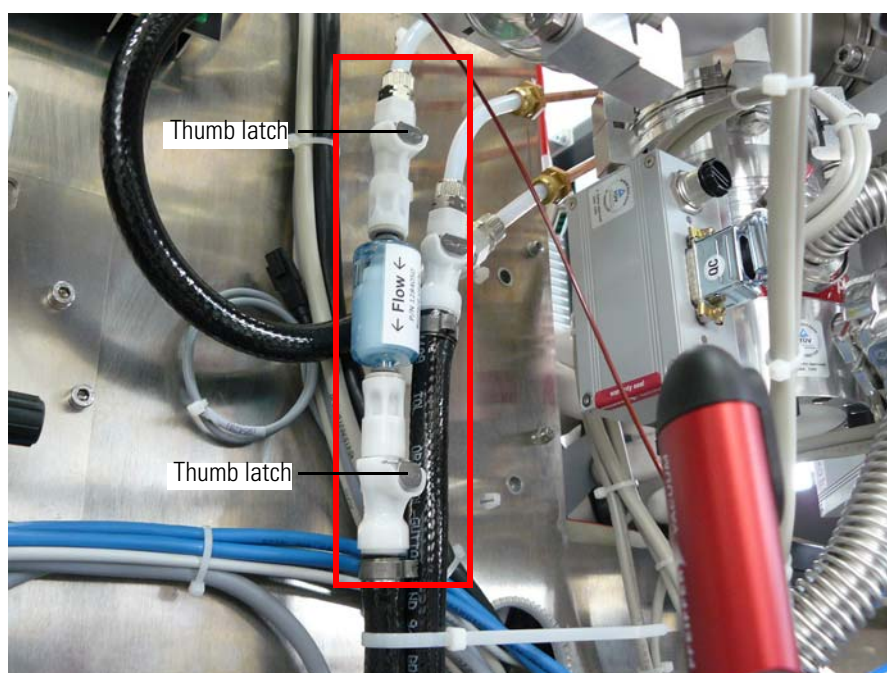


Figure 3-50. Installed water filter

4. Remove the filter from the assembly by using a 5/16 inch wrench and pressing against the gray ring on the quick coupling, away from the filter. See [Figure 3-51](#).

User Maintenance

Maintenance of the Cooling Circuit

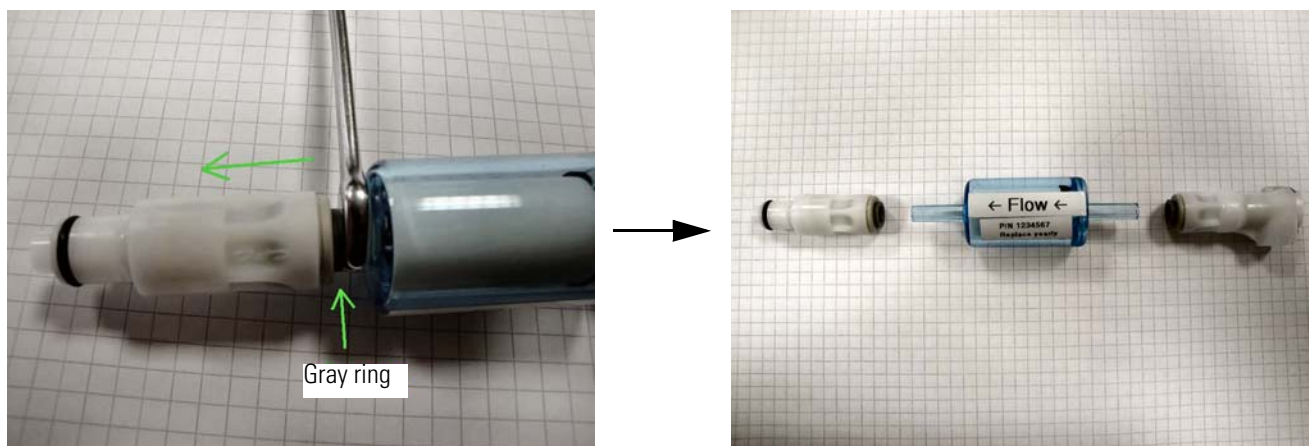


Figure 3-51. Removing the filter cartridge

5. Insert the new filter into the quick couplings. See [Figure 3-52](#).



Figure 3-52. Filter cartridge with Quick couplings

Caution Pay special attention to the direction of flow. Reversing the flow can damage both the flow sensor and the filter. ▲

6. Connect the filter assembly to the hoses of the cooling lines.
7. Switch on the chiller system. Check for leaks and check the water level in the chiller. Refill as appropriate.
8. Set the instrument to operating condition.

Chapter 4 Replaceable Parts

This chapter contains part numbers for replaceable and consumable parts for the mass spectrometer, data system, and kits. To ensure proper results in servicing the Orbitrap Velos Pro system, order only the parts listed or their equivalent.

Note Not all parts are available for purchase separately. Some parts may only be available for purchase as part of a kit or assembly. ▲

For information on how to order parts, see “[Contacting Us](#)” in the front section of this guide.

Ion Sources

ESI probe, for Ion Max source	OPTON-20011
Low flow metal needle for API 2 probes	OPTON-30004
Nanospray II Ion Source	OPTON-20050
Static Nanospray	OPTON-20051
Dynamic Nanospray	OPTON-97017
APCI probe	OPTON-20012
APPI probe	OPTON-20026
HESI-II Probe Kit	OPTON 20037
High-flow needle insert assembly	OPTON-53010
Low-flow needle insert assembly	OPTON-53011

Parts for the Basic System

Orbitrap Analyzer Installation Kit

Tube, 1/8" × 2.1- 1.4301	0261000
Ferrule, stainless steel; R. 1/8"	0520950
Ferrule, stainless steel; V. 1/8"	0522520
Cap nut, stainless steel; 1/8"	0520890
Fitting KJH06-00.	1221620

Pumping System

For a schematical overview of the pumping system, see [Figure 1-22](#) on [page 1-30](#).

T-piece 13 mm	0512360
Hose 13 × 3.5; PVC	0690720
Turbomolecular pump HiPace 300.	1272910
Turbomolecular pump; HiPace 80	1272920
UHV gauge IKR 270; short	1181380
Compact Pirani Gauge TPR280.	1156400
Water cooling for HiPace 300.	1272930
Water cooling for HiPace 80.	0794742
PVC hose, with steel helix; ID=45 mm, L=1.6 m.	1184330
Hose nipple, DN 40, ISO-KF-45.	1159230
Venting flange; DN 10, KF-G1/8"	1184400
Splinter shield for TMPs, with DN 100 CF-F flange	1198590
Centering ring, with integrated splinter shield; DN 63 ISO	1198600
Anti-magnetic cover for IKR 270	1181390
Gasket; NW 100 CF, copper	0552440
Hose, metal; KF16-KF25 - 250mm	1154130
Gasket; copper, NW 35	0550480
Metal tube, KF NW16×250.	0524260
KF Tee piece; NW 16 KF, stainless steel	0524230
Centering ring with o-ring; DN 16, Viton	0522140
Centering ring with o-ring; DN 25, Viton	0522150
Centering ring; NW 16/10, aluminum-Viton	0522200
Clamping ring; NW 10/16, KF	0521830
Clamping ring; NW 20/25, KF	0521560
Reducing cross piece; DN40/DN16 KF.	1184310
Metal tube; DN40x500	1184350
Metal tube; DN40x750	1181290
Hose clamp; NW 40	1181320
Centering ring; NW 40 KF, aluminum-Viton	0522260
Tension ring; NW 32/40 KF	1181250
Clamping screw; DN63-100 ISO, aluminum.	1042670
Flexible metal hose KF NW 16x500.	0534500

Water Supply

For a schematical overview of the cooling water circuit, see [Figure 1-31](#) on [page 1-42](#).

Quick coupling insert; 9.6 mm	1141640
Quick coupling body; 9.6 mm.	1138960
Hose; 9 x 3, black, PVC	1049540
Hose; 6 x 1, Teflon	1042660
Quick coupling insert; Delrin Acetal, NW 6.4	1185030
Quick coupling body; Delrin Acetal, NW 6.4.....	1185020
Clamping piece 8/16	0370130
Adaptor hose nipple; male, 1/2 x 10	1185840
Flow control sensor	1191740
Filter cartridge; 50 µm DIF	1284050

Gas Supply

For a schematical overview of the gas supply, see [Figure 1-27](#) on [page 1-37](#).

Bulkhead union; 1/16", for hose 4 x 1 (for P/N 069 1130)	1153660
Bulkhead union; 1/8"×1/8"	0523450
Hose; 4 x 1, Teflon	0690280
Hose; 4 x 1, polyurethane, blue	0691130
Capillary 1/16" ID-SS	0605470
Plug-in T-piece; 3 x 6mm	1128140
Regulator + manometer f. Orbitrap	1257670
Capillary; 1/16"x0.13x400mm (red), PEEK	1253830
Coupling; 1/16", SS-100-6	0524340
Ferrule; 1/16" GVF 16-000.....	1121110
Reducer Swagelok; 1/8" × 1/16", stainless steel	0662880
Ferrule; 1/16" GVF/16	0674800
Connector 1/8", for hose OD 4 mm	1128680
Cap nut; 1/16", stainless steel	0520880
Hose; 2 x 1, PTFE.....	1091650
Sleeve; Ø 6 mm.....	1047320
Capillary; PEEKsil, 1/16", 0.1 × 500 mm	1223420
Plug, KQ2P-06	1185620
Cap, KQ2C-06	1258220

Parts Lists for the ETD System

This section contains parts lists for the components of the ETD System of the Orbitrap Velos Pro ETD mass spectrometer.

Quadrupole Orbitrap analyzer, complete	1239200
Housing HCD/ETD	1231740
Separating plate HCD/ETD	1231770
Screw M 4 x 8 DIN912	1044420
O-ring 129 X 4 A Viton	1240520
Lid HCD/ETD	1231760
Screw-in connector; 1/16"	1186150
Gasket; NW 63 ISO, aluminum/Viton	0554060
Blank flange; stainless steel, NW 63	0652620
Clamping screw; DN63-100, aluminum	1028380
Washer 8.4; stainless steel	0470070
Screw, hexagonal; M 8 x 35, stainless steel	0454400
Washer 8.4 x 11 x 1.5, stainless steel	0470860
Screw M8 x 35; stainless steel	0454250
Centering ring NW 16 Viton	0522140
Feedthrough; 8-fold 1,5kV DN16KF	1231750
O-ring; 118 X 5 A, Viton	1168240
Box f. feedthrough; KF16 / Sub-D9	1231800
Cylindrical Screw ISO4762-M6X12-A4	0453300

Vacuum system ETD

Flange clamp; KF16	1145860
PUMP, TURBO, EDWARDS EXT75DX ISO100, TNR	00108-01-00016
Splinter guard, DN_63_ISO	1198600
Metal hose, DN 16 ISO-KF x500	1181410
Centering ring with o-ring; DN 16, Viton	0522140
Clamping ring; NW 10/16, KF	0521830
Hose flange; NW 25 KF	1042330
Centering ring, with o ring; DN 25, Viton	0522150
Clamping ring, DN 25	0521560
Flange, KF16 - hose OD 19	1239340
Reducer; DN 16/DN 25, aluminum	0522160
PUMP, ROTARY VANE, EDWARDS RV3	00108-01-0008
KIT, ACCESSORY, MECHANICAL PUMP, RV3	98000-620007

Gas Supply ETD

Plug-in T-piece; 3 x 6mm	1128140
Sleeve 6 mm	1047320
Reducing hose connector, 3.2->6	1239220
Hose cutter	1239280
T-piece 13 mm	0512360
Hose 9 X 3; PVC, black	1049540
Clamping piece 8/16	0370130

Replaceable Parts

Parts Lists for the ETD System

Hose; 4 X 1, Teflon	0690280
Clamping ring; stainless steel, NW10/16.	1149200
PEEK capillary; 1/16" x 0.040.	1245940
Ferrule, 1/16", for GVF/16	0674800
Ferrule 1/16" - CTFE, collapsible	1224700

Electronic Parts ETD

Cable Y-ADAPTER/T.PUMP	2108630
Coupling; RJ45 BU/2BU	2075210
Patch cable; 0.51MT RJ45 gray SFTP	2080870
Cable POWER DIS./T-PUMP	2081200
Cable CLT-OFFSET-A	2108710
Cable IOS ETD/ION OPTIC-S	2108820
UNIT_ION OPTIC SUPPLY ETD	2108920
PCB LTQ CABLE DRIVER.	2097780
PCB ORBITRAP CABLE RECEIVER.	2097830
PCB ETD CABLE RECEIVER	2097800
Cable ETD/LTQ/ORBITRAP-INTERCONNECT 60	2108940
Cable ETD/LTQ/ORBITRAP-INTERCONNECT 36	2108950
Cable ETD/LTQ/ORBITRAP-INTERCON SUPPLY	2108960
Coaxial cable; ETD IOS/ANALOG CTRL, J5554	2108990
Coaxial cable; ETD IOS/ANALOG CTRL, J5555	2109000
Coaxial cable; ETD IOS/ANALOG CTRL, J5556	2109010
Cable ANALOG CTRL / ETD IOS	2108970
Cable ETD IOS/HCD multipole	2108980
Extension cable; 16A C20-C19 2M	2097050
ADAPTER_IOS/MULTIPOLE-HCD	2100410

Reagent Inlet Module

ASSY, TOOL, ION VOLUME INSERTION/REMOVAL . . .	98000-60028
TUBING, PEEKsil, 1/16" OD, 100mm LONG, RoHS. . .	00109-02-00020
FRLE, 1HOLE, 1/16 OD, 0.4mm ID, VESP/GRPHT, RO	00101-08-00005
FRLE, 2HOLS, 1/16 OD, 0.4mm ID, VESP/GRPHT, RO	00101-08-00006

Inlet Valve Seal Kit. 119265-0003

Inlet valve seal removal tool.	119283-0001
Spool inlet valve seal	119683-0100
O-ring Viton, 0530 ID x 0.082 W	3814-6530

ETD Reagent Kit

ETD Reagent Kit	98000-62008
Angiotensin I, 1mg	00301-15517
Fluoranthene, 150mg	00301-01-0013

The fluoranthene in your ETD Reagent Kit is Sigma/Aldrich Supelco #48535. The fluoranthene MSDS is obtained from the MSDS link at:

www.sigmaaldrich.com/catalog/search/ProductDetail/SUPELCO/48535

Thermo Fisher Scientific supplies fluoranthene as a two vial kit. One vial contains 150 mg of fluoranthene and the other is the required empty vial.

The angiotensin I in your ETD Reagent Kit is Angiotensin I human acetate hydrate (Sigma/Aldrich #A9650). Angiotensin I is potentially hazardous. Handle it in accordance with its MSDS. The angiotensin I MSDS is obtained from the MSDS link at:

www.sigmaaldrich.com/catalog/search/ProductDetail/SIGMA/A9650

Appendix A Fluoranthene

Fluoranthene is used as the Electron Transfer Dissociation (ETD) reagent in the ETD Module portion of the Orbitrap Velos Pro ETD mass spectrometer. The fluoranthene radical anion is generated according to the reaction shown in Figure A-1.

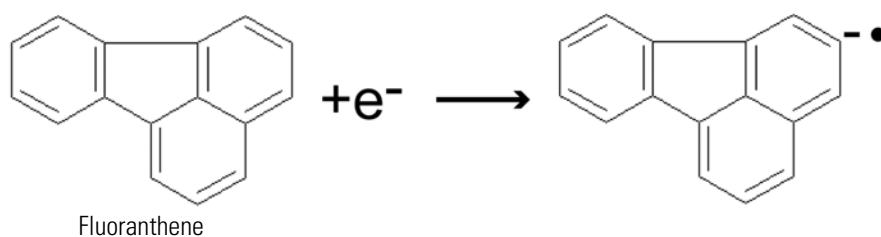


Figure A-1. ETD Reagent (fluoranthene radical anion) generation from fluoranthene

Fluoranthene is potentially hazardous. Use it in accordance with its Material Safety Data Sheet (MSDS).

Note Store and handle all chemicals in accordance with standard safety procedures. The MSDS describing the chemicals being used should be freely available to lab personnel for them to examine at any time. Material Safety Data Sheets (MSDSs) provide summarized information on the hazard and toxicity of specific chemical compounds. The MSDS also provides information on the proper handling of compounds, first aid for accidental exposure, and procedures for cleaning spills or dealing with leaks. Producers and suppliers of chemical compounds are required by law to provide their customers with the most current health and safety information in the form of an MSDS. Read the MSDS for each chemical you use. Dispose of all laboratory reagents in the appropriate way (see the MSDS). ▲

The fluoranthene contained in the ETD Reagent Kit (P/N 98000-62008, see page 4-7) is Sigma/Aldrich Supelco #48535. The fluoranthene MSDS is obtained from the MSDS link at:

www.sigmaaldrich.com/catalog/search/ProductDetail/SUPELCO/48535

Thermo Fisher Scientific supplies fluoranthene as a two vial kit. One vial contains 150 mg of fluoranthene and the other is the required empty vial.

Glossary

This section lists and defines terms used in this guide. It also includes acronyms, metric prefixes, symbols, and abbreviations.

A B C D E F G H I J K L M N O P Q R S T U V W X Y Z

- A**
- A** ampere
- AC** alternating current
- ADC** analog-to-digital converter; a device that converts data from analog to digital form.
- adduct ion** An ion formed by the joining together of two species, usually an ion and a molecule, and often within the ion source, to form an ion containing all the constituent atoms of both species.
- AGC™** See [Automatic Gain Control™ \(AGC\)](#).
- APCI** See [atmospheric pressure chemical ionization \(APCI\)](#).
- APCI corona discharge current** The ion current carried by the charged particles in the APCI source. The voltage on the APCI corona discharge needle supplies the potential required to ionize the particles. The APCI corona discharge current is set; the APCI corona discharge voltage varies, as required, to maintain the set discharge current.
- See also [corona discharge](#) and [APCI corona discharge voltage](#).
- APCI corona discharge voltage** The high voltage that is applied to the corona discharge needle in the APCI source to produce the APCI corona discharge. The corona discharge voltage varies, as required, to maintain the set APCI spray current.
- See also [APCI spray current](#).
- APCI manifold** The manifold that houses the APCI sample tube and nozzle, and contains the plumbing for the sheath and auxiliary gas.
- APCI needle, corona discharge** A needle to which a sufficiently high voltage (typically ± 3 to ± 5 kV) is applied to produce a chemical ionization plasma by the corona discharge mechanism.
- See also [chemical ionization \(CI\)](#), [chemical ionization \(CI\) plasma](#), [atmospheric pressure chemical ionization \(APCI\)](#), and [corona discharge](#).
- APCI nozzle** The nozzle in the APCI probe that sprays the sample solution into a fine mist.
- See also [atmospheric pressure chemical ionization \(APCI\)](#).
- APCI sample tube** A fused silica tube that delivers sample solution to the [APCI nozzle](#). The APCI sample tube extends from the sample inlet to the APCI nozzle.
- See also [atmospheric pressure chemical ionization \(APCI\)](#), and [API stack](#).
- APCI source** Contains the APCI probe assembly, APCI manifold, and API stack.
- See also [atmospheric pressure chemical ionization \(APCI\)](#), [APCI manifold](#), and [API stack](#).
- APCI spray current** The ion current carried by the charged particles in the APCI source. The [APCI corona discharge voltage](#) varies, as required, to maintain the set spray current.
- APCI vaporizer** A heated tube that vaporizes the sample solution as the solution exits the sample tube and enters the atmospheric pressure region of the APCI source.
- See also [atmospheric pressure chemical ionization \(APCI\)](#).

Glossary: API—atmospheric pressure chemical ionization (APCI)

API See [atmospheric pressure ionization \(API\)](#).

API atmospheric pressure region The first of two chambers in the API source. Also referred to as the spray chamber.

API capillary-skimmer region The area between the capillary and the skimmer, which is surrounded by the tube lens. It is also the area of first-stage evacuation in the API source.

API heated capillary A tube assembly that assists in desolvating ions that are produced by the ESI or APCI probe.

See also [API heated capillary voltage](#).

API heated capillary voltage The DC voltage applied to the heated capillary. The voltage is positive for positive ions and negative for negative ions.

See also [API source](#) and [API heated capillary](#).

API ion transfer capillary A tube assembly that assists in desolvating ions that are produced by the ESI, NSI, or APCI probe.

See also [API ion transfer capillary offset voltage](#) and [API ion transfer capillary temperature](#).

API ion transfer capillary offset voltage A DC voltage applied to the ion transfer capillary. The voltage is positive for positive ions and negative for negative ions.

See also [API source](#) and [API ion transfer capillary](#).

API ion transfer capillary temperature The temperature of the ion transfer capillary, which should be adjusted for different flow rates.

See also [API source](#) and [API ion transfer capillary](#).

API source The sample interface between the LC and the mass spectrometer. It consists of the API probe (ESI or APCI) and API stack.

See also [atmospheric pressure ionization \(API\)](#), [ESI source](#), [APCI source](#), [ESI probe](#), and [API stack](#).

API spray chamber The first of two chambers in the API source. In this chamber the sample liquid exits the probe and is sprayed into a fine mist (ESI or NSI) or is vaporized (APCI) as it is transported to the entrance end of the ion transfer capillary.

API spray shield A stainless steel, cylindrical vessel that, in combination with the ESI or APCI flange, forms the atmospheric pressure region of the API source.

See also [atmospheric pressure ionization \(API\)](#).

API stack Consists of the components of the API source that are held under vacuum and includes the [API spray shield](#), [API ion transfer capillary](#), [API tube lens](#), [skimmer](#), the ion transfer capillary mount, and the tube lens and skimmer mount.

See also [atmospheric pressure ionization \(API\)](#) and [API source](#).

API tube lens A lens in the API source that separates ions from neutral particles as they leave the ion transfer capillary. A potential applied to the tube lens focuses the ions toward the opening of the skimmer and helps to dissociate adduct ions.

See also [API tube lens offset voltage](#), [API source](#), [API ion transfer capillary](#), and [adduct ion](#).

API tube lens and skimmer mount A mount that attaches to the heated capillary mount. The tube lens and skimmer attach to the tube lens and skimmer mount.

API tube lens offset voltage A DC voltage applied to the tube lens. The value is normally tuned for a specific compound.

See also [API tube lens](#), [adduct ion](#), and [source CID](#).

AP-MALDI See [atmospheric pressure matrix-assisted laser desorption/ionization \(AP-MALDI\)](#).

APPI See [Atmospheric Pressure Photoionization \(APPI\)](#).

ASCII American Standard Code for Information Interchange

atmospheric pressure chemical ionization (APCI) A soft ionization technique done in an ion source operating at atmospheric pressure. Electrons from a corona discharge initiate the process by ionizing the mobile phase vapor molecules. A reagent gas forms, which efficiently produces positive and negative ions of the analyte through a complex series of chemical reactions.

See also [electrospray ionization \(ESI\)](#).

atmospheric pressure ionization (API) Ionization performed at atmospheric pressure by using [atmospheric pressure chemical ionization \(APCI\)](#), [electrospray ionization \(ESI\)](#), or [nanospray ionization \(NSI\)](#).

atmospheric pressure matrix-assisted laser desorption/ionization (AP-MALDI) Matrix-assisted laser desorption/ionization in which the sample target is at atmospheric pressure.

See also [matrix-assisted laser desorption/ionization \(MALDI\)](#).

Atmospheric Pressure Photoionization (APPI) A soft ionization technique in which an ion is generated from a molecule when it interacts with a photon from a light source.

atomic mass unit Atomic Mass Unit (u) defined by taking the mass of one atom of carbon-12 as being 12u; unit of mass for expressing masses of atoms or molecules.

Automatic Gain Control™ (AGC) Sets the ion injection time to maintain the optimum quantity of ions for each scan. With AGC on, the scan function consists of a prescan and an analytical scan.

See also [ion injection time](#).

auxiliary gas The outer-coaxial gas (nitrogen) that assists the sheath (inner-coaxial) gas in dispersing and/or evaporating sample solution as the sample solution exits the APCI, ESI, or H-ESI nozzle.

auxiliary gas flow rate The relative rate of flow of [auxiliary gas](#) (nitrogen) into the API source reported in arbitrary units.

auxiliary gas inlet An inlet in the API probe where auxiliary gas is introduced into the probe.

See also [auxiliary gas](#) and [atmospheric pressure ionization \(API\)](#).

auxiliary gas plumbing The gas plumbing that delivers outer coaxial nitrogen gas to the ESI or APCI nozzle.

auxiliary gas valve A valve that controls the flow of auxiliary gas into the API source.

B

b bit

B byte (8 b)

baud rate data transmission speed in events per second

BTU British thermal unit, a unit of energy

C

°C degrees Celsius

CE central electrode (of the Orbitrap analyzer);

European conformity. Mandatory European marking for certain product groups to indicate conformity with essential health and safety requirements set out in European Directives.

cfm cubic feet per minute

chemical ionization (CI) The formation of new ionized species when gaseous molecules interact with ions. The process can involve transfer of an electron, proton, or other charged species between the reactants.

chemical ionization (CI) plasma The collection of ions, electrons, and neutral species formed in the ion source during chemical ionization.

See also [chemical ionization \(CI\)](#).

CI See [chemical ionization \(CI\)](#).

CID See [collision-induced dissociation \(CID\)](#).

cm centimeter

cm³ cubic centimeter

collision gas A neutral gas used to undergo collisions with ions.

collision-induced dissociation (CID) An ion/neutral process in which an ion is dissociated as a result of interaction with a neutral target species.

consecutive reaction monitoring (CRM) scan type A scan type with three or more stages of mass analysis and in which a particular multi-step reaction path is monitored.

Convectron™ gauge A thermocouple bridge gauge that is sensitive to the pressure as well as the thermal conductivity of the gas used to measure pressures between X and Y.

corona discharge In the APCI source, an electrical discharge in the region around the corona discharge needle that ionizes gas molecules to form a chemical ionization (CI) plasma, which contains CI reagent ions.

See also [chemical ionization \(CI\) plasma](#) and [atmospheric pressure chemical ionization \(APCI\)](#).

CPU central processing unit (of a computer)

CRM See [consecutive reaction monitoring \(CRM\) scan type](#).

C-Trap curved linear trap

<Ctrl> control key on the terminal keyboard

D

d depth

Da dalton

DAC digital-to-analog converter

damping gas Helium gas introduced into the ion trap mass analyzer that slows the motion of ions entering the mass analyzer so that the ions can be trapped by the RF voltage fields in the mass analyzer.

data-dependent scan A scan mode that uses specified criteria to select one or more ions of interest on which to perform subsequent scans, such as MS/MS or ZoomScan.

DC direct current

divert/inject valve A valve on the mass spectrometer that can be plumbed as a divert valve or as a loop injector.

DS data system

DSP digital signal processor

E

ECD See [electron capture dissociation \(ECD\)](#).

EI electron ionization

electron capture dissociation (ECD) A method of fragmenting gas phase ions for tandem mass spectrometric analysis. ECD involves the direct introduction of low energy electrons to trapped gas phase ions.

See also [electron transfer dissociation \(ETD\)](#) and [infrared multiphoton dissociation \(IRMPD\)](#).

electron multiplier A device used for current amplification through the secondary emission of electrons. Electron multipliers can have a discrete dynode or a continuous dynode.

electron transfer dissociation (ETD) A method of fragmenting peptides and proteins. In electron transfer dissociation (ETD), singly charged reagent anions transfer an electron to multiply protonated peptides within the ion trap mass analyzer. This leads to a rich ladder of sequence ions derived from cleavage at the amide groups along the peptide backbone. Amino acid side chains and important modifications such as phosphorylation are left intact.

See also [fluoranthene](#).

electrospray ionization (ESI) A type of atmospheric pressure ionization that is currently the softest ionization technique available to transform ions in solution into ions in the gas phase.

EMBL European Molecular Biology Laboratory

<Enter> Enter key on the terminal keyboard

ESD ElectroStatic Discharge. Discharge of stored static electricity that can damage electronic equipment and impair electrical circuitry, resulting in complete or intermittent failures.

ESI See [electrospray ionization \(ESI\)](#).

ESI flange A flange that holds the [ESI probe](#) in position next to the entrance of the heated capillary, which is part of the API stack. The ESI flange also seals the atmospheric pressure region of the API source and, when it is in the engaged position against the spray shield, compresses the high-voltage safety-interlock switch.

ESI probe A probe that produces charged aerosol droplets that contain sample ions. The ESI probe is typically operated at liquid flows of 1 $\mu\text{L}/\text{min}$ to 1 mL/min without splitting. The ESI probe includes the ESI manifold, sample tube, nozzle, and needle.

ESI source Contains the ESI probe and the API stack.

See also [electrospray ionization \(ESI\)](#), [ESI probe](#), and [API stack](#).

ESI spray current The flow of charged particles in the ESI source. The voltage on the ESI spray needle supplies the potential required to ionize the particles.

ESI spray voltage The high voltage that is applied to the spray needle in the ESI source to produce the ESI spray current. In ESI, the voltage is applied to the spray liquid as it emerges from the nozzle.

See also [ESI spray current](#).

ETD See [electron transfer dissociation \(ETD\)](#).

eV Electron Volt. The energy gained by an electron that accelerates through a potential difference of one volt.

Extensible Markup Language See [XML \(Extensible Markup Language\)](#).

external lock mass A lock that is analyzed in a separate MS experiment from your sample. If you need to run a large number of samples, or if accurate mass samples will be intermingled with standard samples, you might want to use external lock masses. These allow more rapid data acquisition by eliminating the need to scan lock masses during each scan.

See also [internal lock mass](#).

F

f femto (10^{-15})

°F degrees Fahrenheit

.fasta file extension of a SEQUEST™ search database file

ft foot

Fast Fourier Transform (FFT) An algorithm that performs a Fourier transformation on data. A Fourier transform is the set of mathematical formulae by which a time function is converted into a frequency-domain function and the converse.

FFT See [Fast Fourier Transform \(FFT\)](#).

fluoranthene A reagent anion that is used in an [electron transfer dissociation \(ETD\)](#) experiment.

firmware Software routines stored in read-only memory. Startup routines and low-level input/output instructions are stored in firmware.

forepump The pump that evacuates the foreline. A rotary-vane pump is a type of forepump.

Fourier Transform - Ion Cyclotron Resonance Mass Spectrometry (FT-ICR MS) A technique that determines the mass-to-charge ratio of an ion by measuring its cyclotron frequency in a strong magnetic field.

fragment ion A charged dissociation product of an ionic fragmentation. Such an ion can dissociate further to form other charged molecular or atomic species of successively lower formula weights.

fragmentation The dissociation of a molecule or ion to form fragments, either ionic or neutral. When a molecule or ion interacts with a particle (electron, ion, or neutral species) the molecule or ion absorbs energy and can subsequently fall apart into a series of charged or neutral fragments. The mass spectrum of the fragment ions is unique for the molecule or ion.

FT Fourier Transformation

FT-ICR MS See [Fourier Transform - Ion Cyclotron Resonance Mass Spectrometry \(FT-ICR MS\)](#).

FTMS Fourier Transformation Mass Spectroscopy

full-scan type Provides a full mass spectrum of each analyte or parent ion. With the full-scan type, the mass analyzer is scanned from the first mass to the last mass without interruption. Also known as single-stage full-scan type.

FWHM Full Width at Half Maximum

G

g gram

G Gauss; giga (10^9)

GC gas chromatograph; gas chromatography

GC/MS gas chromatography / mass spectrometer

GUI graphical user interface

H

h hour

h height

handshake A signal that acknowledges that communication can take place.

HCD See [higher energy collision-induced dissociation \(HCD\)](#).

header information Data stored in each data file that summarizes the information contained in the file.

H-ESI probe Heated-electrospray ionization (H-ESI) converts ions in solution into ions in the gas phase by using [electrospray ionization \(ESI\)](#) in combination with heated [auxiliary gas](#).

higher energy collision-induced dissociation (HCD)

Collision-induced dissociation that occurs in the HCD cell of the [Orbitrap mass analyzer](#). The HCD cell consists of a straight multipole mounted inside a collision gas-filled tube. A voltage offset between C-Trap and HCD cell accelerates parent ions into the collision gas inside the HCD cell, which causes the ions to fragment into product ions. The product ions are then returned to the Orbitrap analyzer for mass analysis. HCD produces triple quadrupole-like product ion mass spectra.

high performance liquid chromatography (HPLC)

Liquid chromatography in which the liquid is driven through the column at high pressure. Also known as high pressure liquid chromatography.

HPLC See [high performance liquid chromatography \(HPLC\)](#).

HV high voltage

Hz hertz (cycles per second)

I

ICR ion cyclotron resonance

ID inside diameter

IEC International Electrotechnical Commission

IEEE Institute of Electrical and Electronics Engineers

in. inch

infrared multiphoton dissociation (IRMPD) In infrared multiphoton dissociation (IRMPD), multiply charged ions consecutively absorb photons emitted by a infrared laser until the vibrational excitation is sufficient for their fragmentation. The fragments continue to pick up energy from the laser pulse and fall apart further to ions of lower mass.

See also [electron capture dissociation \(ECD\)](#).

instrument method A set of experiment parameters that define Xcalibur operating settings for the autosampler, liquid chromatograph (LC), mass spectrometer, divert valve, syringe pump, and so on. Instrument methods are saved as file type .meth.

internal lock mass A lock that is analyzed during the same MS experiment as your sample and is contained within the sample solution or infused into the LC flow during the experiment. Internal lock masses provide the most accurate corrections to the data.

See also [external lock mass](#).

I/O input/output

ion gauge Measures the pressure in the mass analyzer region (high vacuum region) of the vacuum manifold.

ion injection time The amount of time that ions are allowed to accumulate in the ion trap mass analyzer when AGC is off. With AGC on, the ion injection time is set automatically (up to the set maximum ion injection time) based on the AGC target value.

See also: [Automatic Gain Control™ \(AGC\)](#).

ion optics Focuses and transmits ions from the API source to the mass analyzer.

ion source A device that converts samples to gas-phase ions.

ion sweep cone A removable cone-shaped metal cover that fits on top of the [API ion transfer capillary](#) and acts as a physical barrier to protect the entrance of the capillary.

ion sweep gas Extra nitrogen gas that flows along the axis of the API ion transfer capillary (between the ion sweep cone and the capillary block) towards the API spray. The sweep gas flow is thus countercurrent to the flow of the ions.

See also [ion sweep gas pressure](#).

ion sweep gas pressure The rate of flow of the sweep gas (nitrogen) into the API source. A measurement of the relative flow rate (in arbitrary units) to provide the required flow of nitrogen gas out from the Ion Sweep cone towards the API spray.

See also [ion sweep gas](#).

IRMPD See [infrared multiphoton dissociation \(IRMPD\)](#).

K

k kilo (10^3 , 1000)

K kilo (2^{10} , 1024)

KEGG Kyoto Encyclopedia of Genes and Genomes

kg kilogram

L

l length

L liter

LAN local area network

lb pound

LC See [liquid chromatography \(LC\)](#).

LC/MS See [liquid chromatography / mass spectrometry \(LC/MS\)](#).

LED light-emitting diode

LHe liquid helium

liquid chromatography (LC) A form of elution chromatography in which a sample partitions between a stationary phase of large surface area and a liquid mobile phase that percolates over the stationary phase.

liquid chromatography / mass spectrometry (LC/MS) An analytical technique in which a high-performance liquid chromatograph (LC) and a mass spectrometer (MS) are combined.

LN2 liquid nitrogen

lock mass A known reference mass in the sample that is used to correct the mass spectral data in an accurate mass experiment and used to perform a real-time secondary mass calibration that corrects the masses of other peaks in a scan. Lock masses with well-defined, symmetrical peaks work best. You can choose to use [internal lock mass](#) or [external lock mass](#).

log file A text file, with a .log file extension, that is used to store lists of information.

M

μ micro (10^{-6})

m meter; milli (10^{-3})

M mega (10^6)

M⁺ molecular ion

MALDI See [matrix-assisted laser desorption/ionization \(MALDI\)](#).

matrix-assisted laser desorption/ionization (MALDI) A method of ionizing proteins where a direct laser beam is used to facilitate vaporization and ionization while a matrix protects the biomolecule from being destroyed by the laser.

MB Megabyte (1 048 576 bytes)

MH⁺ protonated molecular ion

microscan One mass analysis (ion injection and storage or scan-out of ions) followed by ion detection. Microscans are summed, to produce one scan, to improve the signal-to-noise ratio of the mass spectral data. The number of microscans per scan is an important factor in determining the overall scan time.

min minute

mL milliliter

mm millimeter

MRFA A peptide with the amino acid sequence methionine–arginine–phenylalanine–alanine.

MS mass spectrometer; mass spectrometry

MS MS^n power: where $n = 1$

MS scan modes Scan modes in which only one stage of mass analysis is performed. The scan types used with the MS scan modes are [full-scan type](#) and [selected ion monitoring \(SIM\) scan type](#).

MSDS Material Safety Data Sheet

MS/MS Mass spectrometry/mass spectrometry, or tandem mass spectrometry is an analytical technique that involves two stages of mass analysis. In the first stage, ions formed in the ion source are analyzed by an initial analyzer. In the second stage, the mass-selected ions are fragmented and the resultant ionic fragments are mass analyzed.

MS^n scan mode The scan power equal to 1 to 10, where the scan power is the power n in the expression MS^n . MS^n is the most general expression for the scan mode, which can include the following:

- The scan mode corresponding to the one stage of mass analysis in a single-stage full-scan experiment or a selected ion monitoring (SIM) experiment
- The scan mode corresponding to the two stages of mass analysis in a two-stage full-scan experiment or a selected reaction monitoring (SRM) experiment
- The scan mode corresponding to the three to ten stages of mass analysis ($n = 3$ to $n = 10$) in a multi-stage full-scan experiment or a consecutive reaction monitoring (CRM) experiment

See also [MS scan modes](#) and [MS/MS](#).

multipole A symmetrical, parallel array of (usually) four, six, or eight cylindrical rods that acts as an ion transmission device. An RF voltage and DC offset voltage are applied to the rods to create an electrostatic field that efficiently transmits ions along the axis of the multipole rods.

m/z Mass-to-charge ratio. An abbreviation used to denote the quantity formed by dividing the mass of an ion (in u) by the number of charges carried by the ion. For example, for the ion $C_7H_7^{2+}$, $m/z=45.5$.

N

n nano (10^{-9})

nanospray ionization (NSI) A type of electrospray ionization (ESI) that accommodates very low flow rates of sample and solvent on the order of 1 to 20 nL/min (for static nanospray) or 100 to 1000 nL/min (for dynamic nanospray).

NCBI National Center for Biotechnology Information (USA)

NIST National Institute of Standards and Technology (USA)

NMR Normal Mass Range

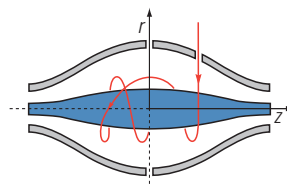
NSI See [nanospray ionization \(NSI\)](#).

O

octapole An octagonal array of cylindrical rods that acts as an ion transmission device. An RF voltage and DC offset voltage applied to the rods create an electrostatic field that transmits the ions along the axis of the octapole rods.

OD outside diameter

Orbitrap mass analyzer The Orbitrap™ mass analyzer consists of a spindle-shaped central electrode surrounded by a pair of bell-shaped outer electrodes. Ions inside the mass analyzer orbit in stable trajectories around the central electrode with harmonic oscillations along it.



Two detection electrodes record an image current of the ions as they undergo harmonic oscillations. A Fourier transformation extracts different harmonic frequencies from the image current. An ion's

mass-to-charge ratio m/z is related to the frequency f of its harmonic oscillations and to the instrumental constant k by:

$$m/z = k/f^2$$

OT Orbitrap

See [Orbitrap mass analyzer](#).

OVC outer vacuum case

Ω ohm

P

p pico (10^{-12})

Pa pascal

parent ion An electrically charged molecular species that can dissociate to form fragments. The fragments can be electrically charged or neutral species. A parent ion can be a molecular ion or an electrically charged fragment of a molecular ion. Also called a precursor ion.

parent mass The mass-to-charge ratio of a parent ion. The location of the center of a target parent-ion peak in mass-to-charge ratio (m/z) units. Also known as precursor mass.

See also: [parent ion](#).

PCB printed circuit board

PDA detector Photodiode Array detector is a linear array of discrete photodiodes on an integrated circuit chip. It is placed at the image plane of a spectrometer to allow a range of wavelengths to be detected simultaneously.

PE protective earth

PID proportional / integral / differential

P/N part number

p-p peak-to-peak voltage

ppm parts per million

PQD pulsed-Q dissociation

precursor ion An electrically charged molecular species that can dissociate to form fragments. The fragments can be electrically charged or neutral species. A precursor ion (PR) can be a molecular ion or an electrically charged fragment of a molecular ion. Also known as parent ion.

precursor mass Mass of the corresponding precursor (or parent) ion or molecule.

psig pounds per square inch, gauge

PTM posttranslational modification

pulsed Q dissociation (PQD) Collision-induced dissociation that involves precursor ion activation at high Q, a time delay to allow the precursor to fragment, and then a rapid pulse to low Q where all fragment ions are trapped. The fragment ions can then be scanned out of the ion trap mass analyzer and detected. PQD eliminates the “1/3 Rule” low mass cut-off for MS/MS data.

Q

quadrupole A symmetrical, parallel array of four hyperbolic rods that acts as a mass analyzer or an ion transmission device. As a mass analyzer, one pair of opposing rods has an oscillating radio frequency (RF) voltage superimposed on a positive direct current (DC) voltage. The other pair has a negative DC voltage and an RF voltage that is 180 degrees out of phase with the first pair of rods. This creates an electrical field (the quadrupole field) that efficiently transmits ions of selected mass-to-charge ratios along the axis of the quadrupole rods.

R

RAM random access memory

raw data Uncorrected liquid chromatograph and mass spectrometer data obtained during an acquisition. Xcalibur and Xcalibur-based software store this data in a file that has a .raw file extension.

resolution The ability to distinguish between two points on the wavelength or mass axis.

retention time (RT) The time after injection at which a compound elutes. The total time that the compound is retained on the chromatograph column.

RF radio frequency

RF lens A multipole rod assembly that is operated with only radio frequency (RF) voltage on the rods. In this type of device, virtually all ions have stable trajectories and pass through the assembly.

RF voltage An AC voltage of constant frequency and variable amplitude that is applied to the ring electrode or endcaps of the mass analyzer or to the rods of a multipole. Because the frequency of this AC voltage is in the radio frequency (RF) range, it is referred to as RF voltage.

RMS root mean square

ROM read-only memory

rotary-vane pump A mechanical vacuum pump that establishes the vacuum necessary for the proper operation of the turbomolecular pump. (Also called a roughing pump or forepump.)

RS-232 An accepted industry standard for serial communication connections. This Recommended Standard (RS) defines the specific lines and signal characteristics used by serial communications controllers to standardize the transmission of serial data between devices.

RT An abbreviated form of the phrase *retention time (RT)*. This shortened form is used to save space when the retention time (in minutes) is displayed in a header, for example, RT: 0.00-3.75.

S

s second

scan mode and scan type combinations A function that coordinates the three processes in the MS detector: ionization, mass analysis, and ion detection. You can combine the various scan modes and scan types to perform a wide variety of experiments.

selected ion monitoring (SIM) scan type A scan type in which the mass spectrometer acquires and records ion current at only one or a few selected mass-to-charge ratio values.

See also [selected reaction monitoring \(SRM\) scan type](#).

selected reaction monitoring (SRM) scan type A scan type with two stages of mass analysis and in which a particular reaction or set of reactions, such as the fragmentation of an ion or the loss of a neutral moiety, is monitored. In SRM a limited number of product ions is monitored.

SEM secondary electron multiplier

Serial Peripheral Interface (SPI) hardware and firmware communications protocol

serial port An input/output location (channel) for serial data transmission.

sheath gas The inner coaxial gas (nitrogen), which is used in the API source to help nebulize the sample solution into a fine mist as the sample solution exits the ESI or APCI nozzle.

sheath gas flow rate The rate of flow of sheath gas into the API source. A measurement of the relative flow rate (in arbitrary units) that needs to be provided at the sheath gas inlet to provide the required flow of [sheath gas](#) to the ESI or APCI nozzle.

sheath gas inlet An inlet in the API probe where [sheath gas](#) is introduced into the probe.

sheath gas plumbing The gas plumbing that delivers [sheath gas](#) to the ESI or APCI nozzle.

sheath gas pressure The rate of flow of sheath gas (nitrogen) into the API source. A measurement of the relative flow rate (in arbitrary units) that needs to be provided at the sheath gas inlet to provide the required flow of inner coaxial nitrogen gas to the ESI or APCI nozzle. A software-controlled proportional valve regulates the flow rate.

See also [sheath gas](#).

sheath gas valve A valve that controls the flow of [sheath gas](#) into the API source. The sheath gas valve is controlled by the data system.

signal-to-noise ratio (S/N) The ratio of the signal height (S) to the noise height (N). The signal height is the baseline corrected peak height. The noise height is the peak-to-peak height of the baseline noise.

SIM See [selected ion monitoring \(SIM\) scan type](#).

skimmer A vacuum baffle between the higher pressure capillary-skimmer region and the lower pressure region. The aperture of the skimmer is offset with respect to the bore of the ion transfer capillary.

source CID A technique for fragmenting ions in an [atmospheric pressure ionization \(API\)](#) source. Collisions occur between the ion and the background gas, which increase the internal energy of the ion and stimulate its dissociation.

SPI See [Serial Peripheral Interface \(SPI\)](#).

SRM See [selected reaction monitoring \(SRM\) scan type](#).

sweep gas Nitrogen gas that flows out from behind the sweep cone in the API source. Sweep gas aids in solvent declustering and adduct reduction.

See also [sweep gas flow rate](#).

sweep gas flow rate The rate of flow of sweep gas into the API source. A measurement of the relative flow rate (in arbitrary units) to provide the required flow of nitrogen gas to the sweep cone of the API source.

See also [sweep gas](#).

syringe pump A device that delivers a solution from a syringe at a specified rate.

T

T Tesla

target compound A compound that you want to identify or quantitate or that a specific protocol (for example, an EPA method) requires that you look for. Target compounds are also called analytes, or target analytes.

TIC See [total ion current \(TIC\)](#).

TMP See [turbomolecular pump](#).

Torr A unit of pressure, equal to 1 mm of mercury and 133.32 Pa.

total ion current (TIC) The sum of the ion current intensities across the scan range in a mass spectrum.

tube lens offset The voltage offset from ground that is applied to the tube lens to focus ions toward the opening of the skimmer.

See also [source CID](#).

Tune Method A defined set of mass spectrometer tune parameters for the ion source and mass analyzer. Tune methods are defined by using the instrument software's tune window and saved as tune file.

A tune method stores tune parameters only. (Calibration parameters are stored separately, not with the tune method.)

tune parameters Instrument parameters whose values vary with the type of experiment.

turbomolecular pump A vacuum pump that provides a high vacuum for the mass spectrometer and detector system.

TWA time weighted average

U

u atomic mass unit

UHV ultra high vacuum

ultra-high performance liquid chromatography (U-HPLC) See [high performance liquid chromatography \(HPLC\)](#).

Ultramark 1621 A mixture of perfluoroalkoxycyclotriphosphazenes used for ion trap calibration and tuning. It provides ESI singly charged peaks at m/z 1022.0, 1122.0, 1222.0, 1322.0, 1422.0, 1522.0, 1622.0, 1722.0, 1822.0, and 1921.9.

UMR Universal Mass Range

V

V volt

VAC volts alternating current

VDC volts direct current

vacuum manifold A thick-walled, aluminum chamber with machined flanges on the front and sides and various electrical feedthroughs and gas inlets that encloses the API stack, ion optics, mass analyzer, and ion detection system.

Glossary: vacuum system—XML (Extensible Markup Language)

vacuum system Components associated with lowering the pressure within the mass spectrometer. A vacuum system includes the vacuum manifold, pumps, pressure gauges, and associated electronics.

vent valve A valve that allows the vacuum manifold to be vented to air or other gases. A solenoid-operated valve.

vol volume

W

w width

W watt

WEEE European Union Waste Electrical and Electronic Equipment Directive. Provides guidelines for disposal of electronic waste.

X

XML (Extensible Markup Language) A general-purpose markup language that is used to facilitate the sharing of data across different information systems, particularly via the Internet.

Index

Numerics

00101-08-00006 3-56
00101-08500 1-38
00101-15500 1-38
00101-2500 1-38
00109-02-00020 3-57
00301-01-0013 4-7
00301-15517 4-7
0690280 1-38
119265-0003 3-45
119283-0001 3-46
119683-0100 3-47
12 pin feedthrough
 harness 3-39
 location 3-34, 3-37
120320-0030 3-42
1275730 3-11
1275740 3-11
1284050 3-58
23827-0008 3-3
23827-0009 3-3
32000-60340 3-14
3814-6530 3-47
98000-20060 3-56
98000-62006 3-56–3-57
98000-62008 4-7, A-1

A

A0301-15101 3-5
Active Pirani gauge (APG) 1-35
active temperature control 1-17
actual vial temperature 3-50
aluminum oxide abrasive, number 600 3-14
analog signals 1-53
analyzer
 chamber temperature 1-62
 system 2-10
 temperature 1-42
angiotensin I 4-7
anti-aliasing filters 1-51
API source
 safety interlock switch 2-10
 settings 2-13
applicators, cotton-tipped 3-14
Automatic Gain Control (AGC) 1-12, 2-15
autosampler 2-8–2-9
auxiliary gas 1-37

axial ion ejection 1-12
axial oscillations
 frequency 1-15–1-16

B

bakeout
 controls 1-6–1-7
 timer 1-7, 3-4–3-5
bakeout procedure
 indication 1-7
 starting 1-6, 3-5
 stopping 1-6, 3-5
base studs 3-41, 3-43
bath gas 1-13
bayonet guide 3-30
bayonet lock 3-30
bayonet pin 3-30
buck/boost transformer 1-10
burnout, of filament 3-54
buttons
 Display Status View 2-11
 On/Standby 2-6
 system bakeout 1-6

C

cable tie 1-9
calibration parameters 2-13
CAN bus 1-55
cartridge heaters 3-41
central electrode
 location 1-14
 power supply box 1-42
 voltages 1-15, 1-65
central electrode power supply board
 cooling 1-17
 diagnostic LEDs 1-66
 location 1-59, 1-65
central electrode pulser board
 diagnostic LEDs 1-62
 location 1-59, 1-61
ceramic lens holder 3-39
ceramic spacer 3-41
charge-sign independent trapping (CSIT) 1-12, 1-21
circuit breakers 1-6–1-7
cleaning
 instrument surface 3-3

- ion source block 3-40
- ion source lens assembly 3-33
- ion volume 3-21
- CLT compartment 1-30
- CLT Offset connector 1-59, 1-64
- CLT RF main board
 - diagnostic LEDs 1-64, 1-66
 - function 1-13
 - location 1-59, 1-63
- CLT RF unit 1-63
- CLT RF voltage 1-13
- Cold Ion Gauge 1-30, 1-32, 1-35–1-36, 1-45, 1-48
- collision cell 1-37
- collision gas
 - C-Trap 1-13
 - HCD 1-13
 - inlet 1-10
 - linear trap 1-38
- Communication LED 1-54
- communication link 1-5
- condenser filter, of recirculating chiller 3-58
- connector, to ETD Interface board 1-22
- control elements
 - inlet valve 1-22
 - instrument 1-5
- control panel 1-6
- control unit, for TMPs 1-32
- Convectron gauge
 - ETD Module 1-22–1-23, 1-25, 1-35
 - linear trap 1-35
- cooling water
 - circuit 1-17, 1-42
 - flow control sensor 1-42
 - properties 1-43
 - temperature 1-42
- cover lid, for bakeout controls 1-7
- C-Trap
 - description 1-13
 - gas supply 1-39
- cycle time 1-4

D

- data acquisition analog board
 - diagnostic LEDs 1-51
 - layout 1-50
 - location 1-49, 1-51
- data acquisition digital PCI board
 - diagnostic LEDs 1-50
 - layout 1-50
 - location 1-49–1-50
- data acquisition unit 1-49
- data communication 1-45
- data system
 - communication 1-5, 2-12

- connection 1-9
- log file 2-3
- DC HV Supply PCB 1-22–1-23, 1-26
- default values, of tune and calibration parameters 2-13
- deflector electrode (DE) voltages 1-65
- Display Status View button 2-11
- distilled water 3-3
- drip pan, of ETD forepump 1-34
- dynamic range, of instrument 1-4

E

- E-clips 3-39
- electric power, for recirculating chiller 1-43
- electrodynamic squeezing 1-15
- electronic boards
 - data acquisition analog 1-51
 - data acquisition digital PCI 1-50
 - instrument control 1-52
 - ion optic supply 1-60
 - left side 1-59
 - power distribution 1-53
 - right side 1-46
- electronic connections, to linear trap 1-45
- emergency shutdown 2-2
- entrance voltage 1-60
- entry housing 3-34
- error messages 1-36, 1-54, 1-56
- ETD
 - main access panel 3-24
 - maintenance 3-12
 - reagent carrier gas supply 1-11
- ETD Control PCB 1-22–1-23, 1-25–1-27
- ETD forepump
 - cabinet 1-22
 - connections 1-23, 1-34
 - electrical cord 1-34
 - power supply 1-24
 - weight 3-10
- ETD Ion Optic Supply board 1-18–1-19, 1-23, 1-46, 1-49
- ETD Module
 - forepump receptacle 1-24
 - Heater Control PCB 1-26
 - Interface board 1-23, 1-25
 - ion gauge 1-23
 - maintenance 3-12
 - power panel 1-24
 - power receptacle 1-24
 - power supply 1-9
 - power switch 1-24
 - service switch 1-24
 - Standby condition 2-4
- ETD power module
 - function 1-23
 - panel 1-23, 1-25

ETD Reagent Kit 4-7
 ETD side access panel
 interlocks 3-55
 removing 3-19
 ETD turbomolecular pump
 connections 1-23, 1-34
 controller 1-35
 function 1-32
 location 1-22
 maintenance 3-11
 Ethernet 1-9–1-10
 evacuating, after a complete shutdown 3-4
 Exactive Installation Kit 1-38
 exhaust
 hose 1-11, 1-34
 system 1-11, 1-33–1-34
 exit voltage 1-60
 external calibration 1-4
 external connections 1-10
 extracting voltage 1-14
 extraction, of ion packets 1-14

F

fan filters 1-22, 3-57
 ferrules, tightening 3-56
 filament
 assembly 3-21, 3-42
 burnout 3-54
 emission current 3-42
 function 1-26, 1-28
 wire 3-44
 filter, for cooling water 1-42
 fingerprints, removing 3-3
 flatapole 1-32
 flow control
 sensor 1-42
 UHP nitrogen 1-23, 1-26
 flow rates
 cooling water 1-42
 HCD collision gas 1-39
 flow restrictors 1-29
 fluid bag filter, of chiller 3-58
 fluoranthene 4-7, A-1
 focusing potentials 1-67
 foreline hose connection 3-34
 forepump electrical cord 1-34
 forepumps
 cabinet 1-2, 1-6, 1-33
 effluent 1-33
 linear trap 1-30–1-31, 1-33
 location 1-6, 1-12, 1-33
 oil level 1-33
 oil mist filters 1-33, 3-11
 power supply 1-33

 switches 1-33
 forevacuum tube 1-34
 front panel 1-3
 front side, LEDs 1-5
 FT Electronics switch
 functions 1-56
 location 1-8
 fused silica tubing
 from lower oven 3-56
 from upper oven 3-56
 in ETD Module 1-29

G

gas ballast 1-33
 gas pressures 2-10
 gas supply
 C-Trap 1-39
 helium 1-10
 nitrogen 1-10
 schematics 1-37
 gas valves 3-56
 gloves 3-2
 ground wires 1-46
 guide bar
 function 1-22, 3-34
 handle 3-23–3-24
 opening 3-23–3-24
 guide bar opening 3-45

H

handling, hot pump oil 3-10
 harmonic oscillations, in Orbitrap 1-15
 hazardous materials 1-33
 HCD
 collision gas port 1-37
 flow rate, of collision gas 1-39
 fragment spectra 1-18
 housing 1-30, 1-32
 HCD collision cell
 description 1-18
 supply voltages 1-46
 heated dual restrictor enclosure 1-29
 heated transfer line 1-29
 heater control 1-56
 Heater Control PCB 1-22–1-23, 1-27
 heater ring 3-41
 heating element 1-42
 helium
 gas capillary 1-39
 gas pressure 1-38
 inlet 1-10, 1-37
 purging the line 2-11

- supply 1-10
- high voltage power supply board
 - diagnostic LEDs 1-68
 - function 1-13
 - layout 1-67
 - location 1-59, 1-66
 - voltages 1-67
- high voltage, at PCBs 1-44

I

- image current 1-15–1-16, 1-48
- inlet port, for UHP nitrogen 1-22
- inlet valve
 - block 3-34
 - knob 3-24, 3-34, 3-45
 - lever 1-22, 3-23–3-24, 3-34, 3-45
 - plug 3-24, 3-34, 3-45
 - seal kit 3-45
 - seal tool 3-46
 - solenoid 1-22, 3-34
- installation kit, for the reagent inlet module 3-56
- instrument
 - controls 1-6
 - diagnostics 1-44
 - forevacuum measurement 1-35
 - high vacuum measurement 1-35
 - operating voltage 1-9
 - parts 4-2
 - rear side 1-6
 - right side 1-9
 - shutdown 1-53
 - switching off 1-9, 2-2
- instrument control board
 - diagnostic LEDs 1-53
 - location 1-46, 1-49, 1-52
 - software status LEDs 1-53
- internal calibration 1-4
- internal computer
 - add-ons 1-50–1-51
 - data communication 1-45
 - location 1-46, 1-49
 - rebooting 2-12
 - timer 1-50
- ion dephasing 1-16
- ion gauge, for ETD Module 1-22
- ion optic supply board
 - diagnostic LEDs 1-60
 - location 1-59–1-60
- ion optics 1-12
- ion oscillation 1-14
- ion packets 1-14
 - shape 1-16
- ion source
 - assembly 3-37
 - block 3-21, 3-39, 3-41, 3-43

- filament 3-41, 3-43
- filament assembly 3-42
- lens assembly 3-39, 3-43
- lenses 3-21
- PCB 3-39, 3-41, 3-43
- schematics 1-23
- springs 3-39
- ion trajectory 1-14
- ion volume
 - alignment arrow 3-30
 - cleaning 3-21
 - key thumbscrew 3-41
 - location 1-23, 1-29
 - temperature 3-28
- ion volume assembly, cleaning frequency 3-21
- ion volume tool
 - alignment arrow 3-30
 - components 3-23
 - handle 1-22, 3-25
 - usage 3-21, 3-27
- ionization techniques 1-12
- ions
 - detection 1-16
 - electrodynamic squeezing 1-15
 - image current 1-16
 - packet shape 1-16–1-17

L

- laboratory
 - exhaust system 1-33–1-34
 - gas supplies 1-10
 - ventilation 1-38
- LC 2-6–2-7
- leak checking 1-26, 1-40
- LEDs
 - power control 1-7
 - system status 1-5, 1-7
 - vacuum 1-5
- left side panel, of instrument 1-3
- lens voltages 1-66
- line power 1-9
- linear trap
 - collision gas 1-38
 - Communication LED 2-9, 2-12
 - connections 1-53
 - description 1-12
 - electronics 1-45
 - forepumps 1-31, 1-33
 - LEDs 1-5
 - location 1-2
 - maintenance 3-1
 - Power LED 2-9, 2-12
 - power panel 1-6, 1-33
 - Reset button 2-12
 - System LED 2-10, 2-12

- turbopump 1-32
- Vacuum LED 2-10
- vacuum measurement 1-35
- vent valve 1-36, 1-39, 2-2–2-3
- Link Port Signal line 1-52
- log file 2-3
- lower oven 3-56

M

- magnet support 3-21
- magnet yoke 3-39
- magnets, in ion source 3-21, 3-39
- main circuit breaker 2-8
- main power switch
 - function 1-9, 1-56
 - location 1-6
 - Off position 2-2
 - securing 1-9
 - status 1-5
- main RF supply 1-63
- mains failure 1-57, 2-2, 3-4
- mains supply
 - for ETD Module 1-9
 - for instrument 1-9
- mains supply, for the ETD Module 1-9
- maintenance
 - API source 1-12
 - cooling circuit 3-58
 - linear trap 3-1
 - procedures 3-2
 - reagent ion source 3-20
 - recirculating chiller 3-58
 - vacuum system 3-4
- mass accuracy 1-4
- mass range 1-4
- Material Safety Data Sheet 3-48
- micro controller 1-52–1-53
- MS/MS 1-4
- MSDS 3-48
- multiple socket outlets 1-7, 1-56, 2-2

N

- nitrogen
 - CLT 1-37
 - connection to source 1-38
 - cooling gas 1-28
 - gas flow 1-39
 - inlet 1-10
 - pressure 1-39
 - pressure regulator 1-39–1-40
 - supply 1-10
 - tubing 1-39
 - venting 1-37

O

- octapole RF voltages 1-60
- oil
 - level 1-33
 - mist filters 1-33–1-34, 3-11
- On/Standby button 2-4, 2-6, 3-22
- operating vacuum 1-5
- Orbitrap
 - applied voltages 1-13
 - central electrode 1-61
 - chamber 1-31
 - control LEDs 1-6
 - detection electrodes 1-61
 - differential pumping 1-14
 - electrodes 1-15
 - Instrument Control board 1-23
 - ion extraction 1-14
 - ion trajectories 1-16
 - layout 1-13
 - lenses 1-14
 - measuring principle 1-14–1-15
 - voltages 1-13, 1-16
- output current, of preamplifier 1-48

P

- padlock, for main power switch 1-9
- parameters
 - calibration 2-13
 - default values 2-13
 - tune 2-13
- part numbers 4-1
- Peltier elements
 - function 1-17, 1-42, 1-62
 - location 1-42, 1-65
 - temperature 1-63
 - water supply 1-32, 1-48
- Penning gauge 1-56
- peripherals power outlet 1-23–1-24
- Pirani gauge 1-30–1-31, 1-36, 1-56
- position, of inlet valve lever 3-34
- power
 - connector 1-6, 1-10
 - connector for linear trap 1-12
 - control LEDs 1-8
 - control panel 1-7–1-8
 - panel 1-6
 - supply for linear trap 1-12
- power distribution
 - operating states 1-56
 - resetting 1-36
 - working modes 1-56
- power distribution board
 - control 1-8
 - function 1-42

- layout 1-54
- location 1-46, 1-53
- power supply 1-7, 1-56
- status LEDs 1-55
- power fail detector 2-3
- Power LED 1-5
- power module 1-22
- power outlet, for peripheral devices 1-9–1-10
- power supply
 - data system 1-9
 - ETD Module 1-9
- power supply 1 board
 - diagnostic LEDs 1-58
 - layout 1-58
 - location 1-46, 1-58
- power supply 2 board
 - diagnostic LEDs 1-52
 - location 1-49
- preamplifier
 - cooling 1-42
 - diagnostic LEDs 1-47–1-48
 - location 1-32, 1-46, 1-48
- pressure readings, in ETD Module 1-35
- pressure regulator
 - location 1-40
 - nitrogen 1-39–1-40
- preventive maintenance 2-8, 3-2
- printed circuit boards 1-44
- proton transfer reactions 1-21
- pumping, the system 2-9, 3-4
- pumps
 - exhaust 1-11
 - forevacuum 1-33
 - manuals 3-5, 3-11
 - oil mist filters 1-33
- purging, helium line 2-11

Q

- quadrupole mass filter 1-21
- quality, of vacuum 1-53

R

- reagent heaters 1-22–1-23, 1-26–1-28
- reagent inlet cover 3-55
- reagent inlet source
 - heating unit 3-51
 - location 3-52
- reagent ion source
 - cleaning frequency 3-13
 - description 1-28–1-29
 - flow restrictors 3-54
 - maintenance 3-20

- maintenance tools 3-14
- Reagent Ion Source dialog box 1-27, 2-4
- Reagent Ion Source instrument control icon 2-4, 3-22
- Reagent Ion Source On check box 1-27–1-28
- reagent vapor, in the carrier gas 1-27
- reagent vials 1-29
- rebooting, of instrument 2-3
- recirculating chiller
 - condenser filter 3-58
 - connections 1-4
 - description 1-43
 - fluid bag filter 3-58
 - maintenance 3-58
- reflector DC voltages 1-60
- removing
 - ETD main access panel 3-18
 - ETD side access panel 3-19
 - gases 2-10, 3-4
 - stains 3-3
- repair covering letter 3-3
- replacing
 - chiller operating fluid 3-58
 - filter cartridge 3-58
 - inlet valve components 3-45
 - ion source filament 3-42
 - operating fluid reservoir of turbopumps 3-11
 - reagent vials 3-48
 - turbopump operating fluid 3-11
- reservoir, of recirculating chiller 3-58
- resetting
 - instrument 2-12
 - system parameters 2-13
 - tune and calibration parameters 2-13
- resolution, of instrument 1-4
- restrictor
 - oven cover 3-55
 - oven heater 1-26
- RF CLT main board 1-66
- RF off & feedback board 1-59, 1-64
- RF output control 1-60
- RF voltage supply 1-63
- right side panel, of instrument 1-3

S

- safety
 - features 1-25
 - interlock switch 2-10
 - problem 3-54
- sample inlet aperture 3-41
- secondary electrons 1-26
- SEM detector 1-12
- serial port connector 1-53
- service switch, of linear trap 1-8
- setting up, conditions for operation 2-10

- sheath gas 1-37
- shutdown 1-57, 2-7, 2-9
- signal communication 1-45
- software
 - debugging 1-53
 - function 1-42, 1-44
 - status LEDs of the instrument control board 1-53
 - Tune Plus window 2-6
- specifications, of instrument 1-4
- spectrum 1-17
- SPI bus
 - function 1-45
 - termination board 1-63–1-64, 1-66, 1-68
- spring clip thumb screw 3-39, 3-41
- stainless steel parts, cleaning 3-14
- stains, removing 3-3
- Standby condition 1-28, 2-4
- starting up
 - ETD Module 2-11
 - instrument 2-9
- status LEDs, of power distribution board 1-55
- Status View 3-22
- sweep gas 1-37
- switches
 - forepumps 1-33
 - FT Electronics 1-8, 3-4
 - linear trap 1-8
 - main power 1-9
 - vacuum pumps 1-8
- switching on, the vacuum system 1-36
- system
 - bakeout 1-5–1-6, 1-36, 1-48, 2-9, 3-4
 - bakeout timer 1-6
 - buttons for bakeout 1-6
 - heating 1-36
 - pump down time 2-10
 - rebooting 2-3
 - resetting 2-12
 - shutdown 2-2
 - Standby 2-4
 - starting up 2-9
 - status LEDs 1-5, 1-54
 - timing 1-50
 - venting 1-39, 2-2
- system status LEDs
 - linear trap 1-2
 - Orbitrap Elite 1-2

T

- temperature
 - analyzer chamber 1-62
 - control 1-17
 - differential 1-17
 - monitoring 3-51

- sensor 3-41
- temperature controller board
 - diagnostic LEDs 1-63
 - function 1-17
 - layout 1-62
 - location 1-59, 1-62
- thermoelectric elements 1-17
- thumbscrews 3-37, 3-39
- tools
 - cleanliness 3-3
 - for cleaning stainless steel parts 3-14
 - for cleaning the ion volume 3-21
 - reagent ion source 3-14
- top lids
 - ETD Module 1-3
 - MS 1-3–1-4
- transfer chamber 1-12
- transfer line
 - bellows 3-37
 - heater 1-26
 - inlet 3-56
 - location 1-23
- transfer multipole 1-19–1-21, 1-32
- triple gas filter 1-41
- tune parameters 2-13
- Tune Plus window
 - buttons 2-6, 2-11
 - diagnostics 1-44
- turbopumps
 - connections 1-53
 - controllers 1-32, 1-35
 - error 1-55
 - HiPace 300 1-30, 1-37, 1-42
 - HiPace 80 1-30, 1-37, 1-42
 - linear trap 1-32, 1-35
 - maintenance 3-11
 - TMP 1 1-30–1-31
 - TMP 2 1-30–1-32
 - TMP 3 1-31–1-32
 - TMP 4 1-31
 - vent valves 1-36

U

- UHP nitrogen
 - inlet port 1-11, 1-40
- UHV chamber
 - components 1-31
 - location 1-30
 - temperature control 1-17
 - vacuum 1-31
- uninterruptible power supply (UPS) 2-3
- upper control panel 1-6–1-7
- user maintenance 3-1–3-2, A-1

V

vacuum

- compartments 1-30
- deterioration 1-36
- failure 1-36, 1-53
- gauges 1-35, 1-53
- quality 1-53
- safety threshold 1-35–1-36
- system 1-9, 1-30

vacuum chamber 1-30–1-31

vacuum components

- left instrument side 1-31
- right instrument side 1-32

Vacuum LED

- instrument 1-5
- linear trap 2-10

vacuum manifold

- location 1-22
- probe plate 3-34–3-35

Vacuum Pumps switch 1-5, 1-8, 1-56

vacuum system

- controls 1-35
- heating 1-36
- moisture 1-36

vent valve

- control 1-39
- function 2-2, 2-7
- linear trap 1-39, 2-3
- nitrogen supply 1-37

ventilation, in the laboratory 1-38

venting pressure 1-40

venting, the system 2-3, 3-19

vial heaters

- cover 3-51–3-52
- location 3-53, 3-56
- ribs 3-53

vial holder 3-52–3-53

vial temperature 1-28

voltage sags 1-10

W

water

- chiller 1-4, 3-58
- filter 1-42, 3-58
- hoses 1-43
- ports 1-10
- temperature 1-43

water cooler, for TMP 1-42

wheels, of instrument 1-3

working modes, of power distribution 1-56

working principle, of the Orbitrap 1-3, 1-20

X

Xcalibur 1-44, 2-15

Thermo Fisher Scientific Inc.

81 Wyman Street

P.O. Box 9046

Waltham, Massachusetts 02454-9046

United States

www.thermoscientific.com