



Nanospray Flex Series Ion Source

User Guide

60053-97127 Revision B • June 2018





© 2018 Thermo Fisher Scientific Inc. All rights reserved.

DirectJunction, EASY-nLC, Exactive, LCQ Deca XP Max, LCQ Fleet, Nanospray Flex, Nanospray Flex NG, nanoViper, and TSQ are trademarks; Unity is a registered service mark; and LCQ, LTQ, Orbitrap, Orbitrap Fusion, Thermo Scientific, Tribrid, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States. Fisher Scientific is a registered trademark of Fisher Scientific Co. in the United States.

The following are registered trademarks in the United States and other countries: Microsoft and Windows are registered trademarks of Microsoft Corporation. Eppendorf is a registered trademark of Eppendorf AG.

The following are registered trademarks in the United States: Dino-Lite and DinoCapture are registered trademarks of AnMo Electronics Corporation. IDEX is a registered trademark of IDEX Corporation. Nanospray is a registered trademark of AB Sciex Pte. Ltd. Swagelok is a registered trademark of Crawford Fitting Company. Tefzel is a registered trademark of The Chemours Company.

SilicaTip is a trademark of New Objective, Inc. GELoader is a trademark of Eppendorf AG.

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

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

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

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

Release history: Rev A, March 2016; Rev B, June 2018

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

WEEE Directive 2012/19/EU



Thermo Fisher Scientific is registered with B2B Compliance (B2Bcompliance.org.uk) in the UK and with the European Recycling Platform (ERP-recycling.org) in all other countries of the European Union and in Norway.

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

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

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

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

Contents

	Preface
	Accessing Documentationix
	Viewing the Product Manuals x
	Viewing Online User Documentation
	Providing Documentation Feedbackx
	Compatible Mass Spectrometers
	Special Notices, Symbols, and Cautionsxi
	Safety Precautions
	Regulatory Compliance
	Contacting Us
Chapter 1	Introduction
•	Advantages of Nanoelectrospray
	Offline Analysis
	Source Housing
Chapter 2	Attaching or Removing the Nanospray Flex Series Ion Source
-	Preparing the Mass Spectrometer
	Attaching the Nanospray Flex Series Ion Source
	Connecting Nitrogen Gas
	Removing the Nanospray Flex Series Ion Source
Chapter 3	Installing the Cameras and LCD Monitor
	Installing the USB Cameras10
	Installing the Analog Cameras and LCD Monitor11
	Adjusting the Video Picture
Chapter 4	Assembling the LC Plumbing13
	Stainless Steel or Glass Emitters
	One- or Two-Column Configuration14
	Using nanoViper Fittings14
	Configuring a Stainless Steel Emitter16
	Configuring a Glass Emitter and Liquid Junction
	Mounting the UHPLC Liquid Junction Cross onto the DirectJunction 21
	Additional Resources

C –

Chapter 5	Configuring the DirectJunction Adapter2	5
-	Attaching the DirectJunction Adapter	5
	Adjusting the Emitter Tip Position	7
Chanter 6	Operating in Offline Nanospray Mode 2	'n
Gliapter o	Office Nersenrey Source Head	.J ^
	Samala Durification	0
	Sample Purification	1
	Attaching the Offline Nanospray Source Head	1 つ
	Installing the Borosilicate Emitter into the Source Head	23
	Opening and Positioning the Borosilicate Emitter	3 4
	Obtaining a Stable Low Flow Spray 3	5
		'
Chapter 7	Configuring the Mass Spectrometer for NSI Mode	7
•	Configuring the NSI Parameters for the Nanospray Flex NG Source	7
	Configuring the NSI Parameters for the Nanospray Flex Source	8
	Selecting the Source for the LCQ Deca XP Max Mass Spectrometer	8
Chapter 8	Troubleshooting4	1
Chanter 9	Renlaceable Parts A	Б
Gliapter 5		5
	Spare rans	s c
	Consumables	0
Appendix A	Contents of the Installation Kits4	7
	Index 4	0
	muex	J

Figures

Figure 1.	Nanospray Flex ion source mounted on a Thermo Scientific mass spectrometer 1
Figure 2.	Nanospray Flex NG ion source (ES072) with the USB cameras (ES218) 3
Figure 3.	Nanospray Flex ion source (ES071) with the USB cameras (ES218)
Figure 4.	Examples of the MS ion sweep cones
Figure 5.	Ion source locking levers (top view) 7
Figure 6.	Nanospray Flex NG ion source connection
Figure 7.	Gas port on the bottom of the ion source
Figure 8.	USB cameras installed on the ion source 10
Figure 9.	Hex screw to secure the top camera11
Figure 10.	nanoViper fitting
Figure 11.	UHPLC fused-silica union (cross section, P/N ES272)21
Figure 12.	Protective cover (side view) 22
Figure 13.	UHPLC liquid junction cross with the bottom protective cover 22
Figure 14.	DirectJunction adapter mounted on the XYZ-manipulator arm (Nanospray
-	Flex)
Figure 15.	Diagram showing the 20-degree emitter angle
Figure 16.	DirectJunction HV connection (left side view) 27
Figure 17.	Adjustment knobs on the XYZ manipulator (front view) 28
Figure 18.	Offline nanospray ion source head (ES260)
Figure 19.	Inserting the StageTip into the borosilicate emitter
Figure 20.	Emitter tip bending on the MS spray cone

F

Figures

Preface

The *Nanospray Flex Series Ion Source User Guide* describes the hardware components, and provides installation and configuration procedures for the Thermo Scientific[™] Nanospray Flex[™] Series ion sources. Depending on the installed camera type, you view the ion source spray by using the camera software or an LCD monitor that is placed on the mass spectrometer (MS).

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

Contents

- Accessing Documentation
- Providing Documentation Feedback
- Compatible Mass Spectrometers
- Special Notices, Symbols, and Cautions
- Safety Precautions
- Regulatory Compliance
- Contacting Us

Accessing Documentation

Thermo Scientific MSs include complete documentation.

- Viewing the Product Manuals
- Viewing Online User Documentation

For system requirements, refer to the release notes on the software DVD.

Viewing the Product Manuals

The Thermo Fisher Scientific service engineer installs the instrument control applications and the instrument manuals on the data system computer.

✤ To view the product manuals

From the Microsoft[™] Windows[™] taskbar, choose **Start > All Apps** (Windows 10) or **All Programs** (Windows 7) **> Thermo Instruments** (or **Thermo** *model*) and so on, where *model* is the specific MS.

Viewing Online User Documentation

Visit the Thermo Fisher Scientific website for product manuals and more.

* To view user documentation from the Thermo Fisher Scientific website

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

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

Providing Documentation Feedback

✤ To suggest changes to the documentation

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



Compatible Mass Spectrometers

Use the Nanospray Flex Series ion source with the appropriate Thermo Scientific MS for nanoelectrospray (commonly referred to as nanoES or nanospray[™]) analysis. Table 1 lists several compatible MSs. For information about your specific instrument, refer to its manuals.



CAUTION If you connect the Nanospray Flex Series ion source to another type of MS, you might impair the safety protection provided by the equipment.

lon source model	Thermo Scientific mass spectrometer
Nanospray Flex NG™	Orbitrap [™] Tribrid [™] Series ^a and TSQ [™] Series II
Nanospray Flex	Exactive [™] Series, LCQ Fleet [™] , LTQ [™] Series, Orbitrap Series, and TSQ Series I

 Table 1.
 Ion sources and compatible mass spectrometers

^a In 2018, the name "Orbitrap Fusion[™] Series" changed to "Orbitrap Tribrid Series."

Special Notices, Symbols, and Cautions

Make sure you understand the special notices, symbols, and caution labels in this guide. Most of the special notices and cautions appear in boxes; those pertaining to safety also have corresponding symbols. Some symbols are also marked on the ion source itself and can appear in color or in black and white. For complete definitions, see Table 2.

Notice, symbol, or label	Meaning
IMPORTANT	Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the product.
Note	Highlights information of general interest.
Тір	Highlights helpful information that can make a task easier.
	Caution: Read the cautionary information associated with this task.

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

Notice, symbol, or label	Meaning
	Chemical hazard: Observe safe laboratory practices and procedures when handling chemicals. Only work with volatile chemicals under a fume or exhaust hood. Wear gloves and other protective equipment, as appropriate, when handling toxic, carcinogenic, mutagenic, corrosive, or irritant chemicals. Use approved containers and proper procedures to dispose of waste oil and when handling wetted parts of the instrument.
	Hot surface: Allow any heated components to cool before touching them.
4	Risk of electric shock: This source uses voltages that can cause electric shock and personal injury. Before removing the source for servicing, shut down the MS and disconnect it from line power.
	Risk of eye injury: Eye injury can occur from splattered chemicals, airborne particles, or sharp objects. Wear safety glasses when handling chemicals or servicing the source.
	Sharp object: Avoid handling the tip of the emitter.
A	Trip obstacle: Be aware of cords or other objects located on the floor.

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

Safety Precautions

Observe the following safety precautions when you operate or perform service on the ion source.



CAUTION Do not perform any servicing other than that contained in this manual. To avoid personal injury or damage to the ion source, do not perform any servicing other than that contained in this manual or related manuals unless you are qualified to do so.



CAUTION The ion source must connect to a certified Thermo Scientific MS, which supplies high voltage (HV) capable of delivering a maximum of 8 kV and 100 μ A. If you connect the source to another type of MS, you might impair the protection provided by the equipment.



CAUTION Be aware of high voltage components. Before you touch a stainless steel or borosilicate emitter, the liquid flowing through a glass emitter, or the offline source head, depending on the plumbing configuration, make sure that you place the MS in off mode.



CAUTION Avoid personal injury. Before you remove the emitter, make sure that you depressurize the LC system. Otherwise, the emitter might eject at a high speed and cause personal injury to you or someone nearby.

Regulatory Compliance

Thermo Fisher Scientific performs complete testing and evaluation of its products to ensure full compliance with applicable North American and European regulations. The Nanospray Flex Series ion sources were tested as part of the Thermo Scientific MS system, which meets the applicable requirements for electromagnetic compatibility (EMC) and product safety. For additional regulatory information, refer to the MS manuals.

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

Ensure continued compliance with regulatory standards:

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

Contacting Us

Contact	Email	Telephone	QR Code ^a					
U.S. Technical Support	us.techsupport.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752						
U.S. Customer Service and Sales	us.customer-support.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752						
Global Support	 To find global contact information or custon 	nize your request	INVER					
	1. Go to thermofisher.com.							
	2. Click Contact Us , select the country, and the you need.	n select the type of support						
	3. At the prompt, type the product name.							
	4. Use the phone number or complete the online form.							
	 To find product support, knowledge bases, and resources 							
	Go to thermofisher.com/us/en/home/technica	d-resources.						
	 To find product information 							
	Go to thermofisher.com/us/en/home/brands/	thermo-scientific.						
Note To provide feedb	back for this document, go to surveymonkey.com/s/	PQM6P62 or send an email	message to					

^a You can use your smartphone to scan a QR Code, which opens your email application or browser.

Technical Publications (techpubs-lcms@thermofisher.com).

1

Introduction

The Nanospray Flex Series ion source, shown in Figure 1, maintains excellent spray stability to ensure efficient evaporation and ionization of the liquid samples—the key to achieving the highest sensitivity at nano-flow rates.

Key benefits of the Nanospray Flex Series are as follows:

- Flexible, user-friendly design (for configuring custom columns)
- Single setup for all online and offline nanoflow applications
- Ability to interface with online nanoscale LC separation techniques

Contents

- Advantages of Nanoelectrospray
- Offline Analysis
- Source Housing



Figure 1. Nanospray Flex ion source mounted on a Thermo Scientific mass spectrometer

Advantages of Nanoelectrospray

The use of electrospray ionization (ESI) has evolved as a leading technique for generating intact, gas-phase ions from thermally labile, polar analytes in solution. In this technique, an emitter (a capillary tube or needle) induces ionization at a controlled distance from a counter electrode. Direct current (dc) voltage is applied, either to the needle or to the solvent, to produce a strong electrical field at the emitter tip. The electric field excites the ions in the solution as they leave the emitter tip. This interaction results in electrohydrodynamic disintegration of the fluid, generation of droplets, and formation of an aerosol jet.

Conventional ESI employs flow rates from 1 μ L/min to 1 mL/min. Expediting desolvation and droplet shrinkage often requires a drying gas, thermal heating, or both, due to the high volume of liquid that exits the emitter. Nanospray ionization (NSI), also known as nanoelectrospray ionization (nanoESI or NSI), is a form of ESI that employs online low rates of 10 to 1000 nL/min. NSI (or nanoESI) generally does not require a drying gas or thermal heating. Compared with ESI, NSI tolerates a wider range of liquid compositions including pure water.

As you lower the flow rate, a lower volume of mobile phase passes through the emitter, producing smaller aerosol droplets. This makes NSI more effective than conventional ESI at concentrating the analyte at the emitter tip, which produces significant increases in sensitivity as demonstrated by the signal response of the MS.

Note The MS's instrument control applications use the terms *nanospray* and *NSI*.

Offline Analysis

Through offline analysis with borosilicate emitters, you can extract the maximum amount of information from very limited amounts of sample. You can also average data extensively to improve the signal-to-noise ratio and conditions optimized for MS/MS experiments. Using the optional offline Nano ES ion source head (ES260), you can do the following:

- Work at low flow rates of 10-40 nL/min.
- Use nearly 100 percent of sample.
- Work effectively with sample volumes down to 300 nL.
- Work with stable sprays over a wide pH range and in high-buffer concentrations.
- Spray from purely aqueous and purely organic solvents.
- Avoid cross-contamination by using disposable emitters.

For additional information, see Operating in Offline Nanospray Mode.

Source Housing

The Nanospray Flex NG (Figure 2) or the Nanospray Flex (Figure 3) ion source is easy to install on the appropriate MS (see Table 1). For installation instructions, see Chapter 2.

The source housing includes two locking levers, an observation cylinder, a position-adjustable column holder, and sliding rails for emitter retraction. Using the observation cylinder, you can view the emitter tip while you move it into position. To enhance your view of the emitter, follow the instructions in Installing the Cameras and LCD Monitor.



Figure 2. Nanospray Flex NG ion source (ES072) with the USB cameras (ES218)



Figure 3. Nanospray Flex ion source (ES071) with the USB cameras (ES218)

2

Attaching or Removing the Nanospray Flex Series Ion Source

Follow these procedures to install or remove the Nanospray Flex Series ion source.

Contents

- Preparing the Mass Spectrometer
- Attaching the Nanospray Flex Series Ion Source
- Connecting Nitrogen Gas
- · Removing the Nanospray Flex Series Ion Source



CAUTION Before you touch a previously operating source or its components, allow the system to cool for a minimum of 20 minutes.

Preparing the Mass Spectrometer

Before you attach the ion source to the instrument, follow this procedure.

To prepare the MS

1. If any other source is installed, remove it after it has cooled to room temperature.

For instructions, refer to the instrument documentation.

2. If the ion sweep cone (Figure 4) is installed, let it cool to room temperature, and then remove it by grasping its outer ridges and pulling it off.

You might need to use a small slotted screwdriver to loosen the screws. The spray cone is directly behind the ion sweep cone (Figure 6).



CAUTION

- Make sure that you do not accidentally lift the release lever located at the top of the API source interface, which could interfere with or prevent installation of the ion source.
- To avoid contaminating the ion transfer tube, do not touch its exposed entrance.





Attaching the Nanospray Flex Series Ion Source

✤ To attach the ion source on the MS

1. Unlock the source's locking levers (Figure 5).



2. Align the two guide pin holes on the back of the source with the guide pins on the front of the MS (Figure 6).





3. Carefully press the source onto the MS, and then lock the locking levers.

Connecting Nitrogen Gas

If you want to provide a controlled environment in the spray area, you can connect nitrogen gas to the ion source.

♦ (Optional) To connect nitrogen gas to the ion source

Connect clean, dry nitrogen gas with a flow rate of 0-5 L/min to the 1/8 in. OD SwagelokTM fitting on the bottom of the source.



Figure 7. Gas port on the bottom of the ion source

HV connector for the DirectJunction™ adapter Swagelok fitting for the gas connection

Removing the Nanospray Flex Series Ion Source

To remove the ion source from the MS

- 1. Turn off the LC liquid flow to the source.
- 2. Disconnect the LC plumbing.
- 3. As applicable, disconnect the cameras and monitor connections, including their power supplies.
- 4. After the system has cooled to room temperature, unlock the source's locking levers.
- 5. Grasp the source housing with both hands and slowly pull it away from the MS.

The MS automatically switches to off mode after a few seconds.

You can now store the source until its next usage. The Nanospray Flex Series ion source does not require cleaning.

3

Installing the Cameras and LCD Monitor

Depending on your camera model (Table 3), follow the applicable procedure to install the two cameras, with or without the monitor. Afterwards, follow the applicable procedure in Adjusting the Video Picture.

Contents

- Installing the USB Cameras
- Installing the Analog Cameras and LCD Monitor
- Adjusting the Video Picture

Table 3. Camera types and installation procedures

Camera	P/N	Monitor?	Procedure ^a
Dino-Lite [™] , USB (2 provided)	ES218	No	To install the USB cameras
Dino-Lite, analog (2 provided)	ES216	Yes ^b	To install the analog cameras and LCD monitor

^a Might require a 1.5 mm hex key to complete.

^b LCD monitor (P/N ES217)

Installing the USB Cameras

Using the DinoCapture[™] application through the data system computer, you can use the monitor to view the spray exiting the emitter tip.

✤ To install the USB cameras

1. Install the camera's DinoCapture application onto the data system computer.

IMPORTANT You must install this application before connecting the camera cables to the data system computer.

2. With the focusing wheel facing toward you, slowly press the cameras into the top and side camera opening (Figure 8).



Figure 8. USB cameras installed on the ion source

- 3. (Optional) Connect each camera to a USB extension cord.
- 4. Plug the USB cords into the data system computer's USB ports.

IMPORTANT Do not connect the cameras to a USB hub.

Tip For troubleshooting purposes, record the USB port used for each camera.

Installing the Analog Cameras and LCD Monitor



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

* To install the analog cameras and LCD monitor

- 1. Install the monitor as follows:
 - a. On the back, attach BNC adapters (if provided) to both Video IN ports.
 - b. Connect the power supply for the monitor between the DC 12V IN socket and an electrical outlet.
 - c. Place the monitor on top of the MS near the source, making sure that it is not too close to the edge of the instrument.
- 2. Install each camera as follows:
 - a. Using a 1.5 mm hex key, loosen the hex screw near the camera opening, and then insert the camera with its focusing wheel facing toward you (Figure 9).
 - b. Tighten the hex screw until it touches the camera-do not overtighten it.







Camera focusing wheel

- c. Connect the yellow video connector to the monitor's Video IN port (see step 1a).
- d. Connect the black connector to its power supply, and then connect the power supply to an electrical outlet.

Adjusting the Video Picture

Depending on your camera model, follow the applicable step to focus the video picture. As needed, refer to the manuals for the cameras and monitor.

To adjust the video picture

- For the USB camera application, do the following:
 - i. In the DinoCapture window, use the mouse to move the visible video image off of the second image, and then adjust the display size as needed.

Tip To learn more about this application, such as the measurement and time-lapse video features, choose **Help > FAQ** or **User's Guide**.

- ii. If the picture is too dark, turn on each camera's light.
- iii. If the picture is out of focus, adjust each camera's focusing wheel.
- For the LCD monitor, do the following:
 - i. Turn on the monitor by pressing its POWER button.
 - ii. If the picture is too dark, turn on each camera's light.
 - iii. If the picture is out of focus, adjust each camera's focusing wheel.
 - iv. To display the other camera input, press the SOURCE button.

The monitor displays the selected video input channel (AV1 or AV2) for a few seconds.

4

Assembling the LC Plumbing

After you decide how many columns you need for your LC analysis and what type of emitter to use, follow the applicable procedure in this chapter to assemble the plumbing for the DirectJunction adapter.



CAUTION Wear protective gloves and safety glasses when handling the solvent lines. To prevent contamination, wear lint-free and powder-free gloves.

Contents

- Stainless Steel or Glass Emitters
- One- or Two-Column Configuration
- Using nanoViper Fittings
- Configuring a Stainless Steel Emitter
- Configuring a Glass Emitter and Liquid Junction
- Mounting the UHPLC Liquid Junction Cross onto the DirectJunction
- Additional Resources

Stainless Steel or Glass Emitters

The MS energizes the stainless steel (SS) emitter through the high-voltage (HV) connection on the ion source's DirectJunction adapter. When you use a glass emitter, the plumbing configuration includes the Liquid Junction adapter, which energizes the liquid flow that passes the HV electrode.

SS emitters are more robust than glass emitters and help maintain a stable spray for longer periods of time.

Glass emitters can have very small openings. However, the very small opening at the emitter tip often results in stability problems associated with blockage, which means that glass emitters rarely last as long as the SS emitters.

Some consider glass emitters to be more bio-inert than SS emitters. This might reduce the risk of nonspecific adsorption of biomolecules and lead to slightly improved sensitivity. However, the degree of this adsorption, and thus the performance change, depends on the chemical characteristics of the sample.

IMPORTANT For glass emitters, do the following:

- Place the Liquid Junction adapter on the high-pressure side of the column. Electrochemical processes that occur at the electrode can otherwise create gas that leads to spray instability.
- Use glass emitters with small ID emitter tips (less than 20 μ m) to create back pressure in the emitter and avoid outgassing and consequent spray instability.

One- or Two-Column Configuration

The Nanospray Flex Series installation kit includes the DirectJunction adapter, which supports both one- and two-column configurations. The following table describes some of the advantages for each configuration.

Configuration	Advantages
One-column	• Fewer connections, which minimize any potential peak broadening that results from dead volumes.
	• Enables MS analysis of compounds that elute during sample loading (as they might not bind to the column material) because the fluid path leads directly to the instrument.
Two-column	• Use of a shorter pre-analytical column (precolumn). This provides an increased loading capacity and loading flow rate when compared to loading sample directly onto the longer analytical column.
	• The precolumn acts as a guard column by protecting the analytical column from particulate matter.

Using nanoViper Fittings

For some LC instruments, the plumbing connections include several nanoViper[™] fittings, such as the one shown in Figure 10. Although these fittings can withstand ultra-high-performance LC (UHPLC) backpressures up to ~1200 bar (~17400 psi), they are fingertight fittings that require only very small torques to seal. Therefore, you must follow the next procedure to avoid damage by overtightening.

Tip Refer to the packing for your nanoViper fittings to determine compatibility with either the 1000 bar capillaries (beige tubing) or the 1200 bar capillaries (blue tubing).

Figure 10. nanoViper fitting



To use a nanoViper fitting

- 1. Insert the nanoViper fitting into the target port and slowly rotate the screw clockwise until you feel resistance.
- 2. Using the black knurled fitting tool, tighten the screw clockwise to an angle of 0–45 degrees (1/8-turn).
- 3. Start operating the system at the desired working pressure and check the backpressure.
- 4. If the backpressure is too low, check the system for leaks.

For instructions, refer to the LC instrument's documentation.

5. If the backpressure continues to be too low, return the system to atmospheric pressure.

IMPORTANT To extend the lifetime of the nanoViper fittings, open and close the connections at only atmospheric system pressures. Opening and closing connections at high system pressures can reduce the lifetime of the fittings.

6. Tighten the screw by as much as an additional 45 degrees. Do not turn the screw beyond an angle of 90 degrees from where you felt the initial resistance.

IMPORTANT To prevent damage to the sealing surface of the nanoViper fitting, do not overtighten the fitting.

Configuring a Stainless Steel Emitter

For a stainless steel (SS) emitter with two columns, follow this HPLC procedure, which shows the recommended configuration. For a one-column configuration, start with Step 2. Make sure that the tubing, sleeves, and unions are compatible and appropriate for the connection method and pressure range.

IMPORTANT Install the emitter last to avoid damaging its tip.

* To configure a stainless steel emitter and two columns for HPLC (max. 300 bar)



To configure a stainless steel emitter and two columns for UHPLC (over 300 bar)



Configuring a Glass Emitter and Liquid Junction

For a glass emitter with two columns, follow the applicable HPLC or UHPLC procedure, which shows the recommended configurations. For a one-column configuration, start with Step 2. Make sure that the tubing, sleeves, and unions are compatible and appropriate for the connection method and pressure range.

IMPORTANT Install the emitter last to avoid damaging its tip.

Note

- Use small ID emitter tips (less than 20 $\mu m)$ to prevent outgassing in the tip and subsequent spray instability.
- The packed-glass emitter contains the stationary phase and has dual roles: as the analytical column and the emitter. The packed-glass emitter is inserted through the ZDV union, which only holds the emitter in place.
- To use a longer analytical column (greater than 20 cm), coil it between the venting Tee and ZDV union.



CAUTION Avoid contact with high voltage. Before you start, make sure that the cable from the UHPLC liquid junction cross is disconnected.

To configure a glass emitter and two columns for HPLC (max. 300 bar)



To configure a glass emitter and two columns for UHPLC (over 300 bar)



CAUTION Avoid contact with high voltage. The UHPLC liquid junction cross operates at high voltage. Make sure that you securely mount the cross to the DirectJunction adapter inside its two-piece protective cover.

IMPORTANT If the columns are less than 360 µm OD, add PEEK sleeves to the ends before inserting them into the UHPLC liquid junction cross.



Mounting the UHPLC Liquid Junction Cross onto the DirectJunction

The UHPLC liquid junction cross, which is included in the UHPLC Liquid Junction Kit (ES269), is for use with glass-emitter LC configurations. Using the tools and equipment listed in Table 4, follow the procedures in this section to mount the cross onto the DirectJunction adapter.

Item	Part number
Gloves, lint-free and powder-free	Fisher Scientific [™] 19-120-2947 ^a
	Unity Lab Services: • 23827-0008 (medium size) • 23827-0009 (large size)
Screwdriver, slotted, small	-
UHPLC Liquid Junction Kit, includes a tightening tool	ES269

Table 4. Tools and equipment

^a Multiple sizes are available.

Figure 11 shows that the ends of the externally threaded PEEK holder have different internal depths. The internal stainless steel cartridge sits within the deeper end of the PEEK holder. The microferrule ends are for use with sleeveless, 360 μ m OD fused-silica tubing, which are secured by the two internally threaded knurled nuts.

Figure 11. UHPLC fused-silica union (cross section, P/N ES272)



* To mount the UHPLC liquid junction cross onto the DirectJunction

1. Using the slotted screwdriver, remove the screws from the top protective cover.

Figure 12 shows a side view of the two-piece protective cover.

Figure 12. Protective cover (side view)



- 2. Place the bottom cover under the DirectJunction's metal rod. If necessary, loosen the compression screw to adjust the hole size, and then tighten the screw.
- 3. Place the UHPLC liquid junction cross into the bottom protective cover, and then attach the top cover by using the slotted screwdriver.

Figure 13. UHPLC liquid junction cross with the bottom protective cover



4. After you complete the plumbing connections, connect the HV cable to the source bottom.

Additional Resources

For general information about configuring Thermo Scientific nanoLC sources, log in (free) to planetorbitrap.com, choose Library > Scientific Library > Keyword, and search for A1969.

Refer to the Thermo Scientific EASY-nLC[™] Series documentation for the following:

- Setting up the column assembly
- Modifying the LC instrument before connecting it to the UHPLC Liquid Junction adapter on the Nanospray Flex Series ion source

For other LC instruments, refer to their documentation.

4 Assembling the LC Plumbing

Additional Resources

Configuring the DirectJunction Adapter

After you attach the DirectJunction adapter to the Nanospray Flex Series ion source, adjust the emitter position.

Contents

- Attaching the DirectJunction Adapter
- Adjusting the Emitter Tip Position

Attaching the DirectJunction Adapter

You can attach the LC plumbing to the DirectJunction adapter before or after you attach the adapter to the source's manipulation arm.

* To attach the DirectJunction adapter to the ion source

- 1. Pull out the bottom of the source to fully expose the XYZ-manipulator arm (Figure 14).
- 2. Using the 3 mm hex key, loosely attach the DirectJunction adapter to the XYZ-manipulator arm with the provided screw.

Note The longer screw provided with the DirectJunction is for the Nanospray Flex NG source, which has a wider manipulator arm.



Figure 14. DirectJunction adapter mounted on the XYZ-manipulator arm (Nanospray Flex) DirectJunction adapter

Bottom of the Hex screw on the XYZ-manipulator arm source

3. Attach the one- or two-column LC configuration to the adapter as shown on page 16, page 19, or page 20.

IMPORTANT Avoid handling the emitter tip when you attach the LC plumbing.

4. Elevate the adapter's free end approximately 20 degrees (Figure 15), and then firmly tighten the screw.

Figure 15. Diagram showing the 20-degree emitter angle





5. Attach the HV cable to the socket on the source's bottom (Figure 16).

The DirectJunction cable (with an SS emitter) or the Liquid Junction cable (with a glass or packed glass emitter) connects to the bottom of the source.

DirectJunction adapter Source bottom HV connector

Figure 16. DirectJunction HV connection (left side view)

Adjusting the Emitter Tip Position

After you attach the DirectJunction Adapter to the manipulator arm, follow this procedure to make any needed positional adjustments for the emitter tip.

✤ To adjust the emitter tip position

- 1. Using the XYZ-manipulator knobs (Figure 17), position the emitter tip almost on-axis with the ion transfer tube.
- 2. As you make adjustments, observe the distance on the monitor. If necessary, follow the procedure, Adjusting the Video Picture.



Figure 17. Adjustment knobs on the XYZ manipulator (front view)

Fine-adjustment for the Zaxis



CAUTION Be aware of accessible high voltage components. Although the source is shielded, you can easily access its head. Always make sure that the MS is in off mode before you touch the source head. Do not leave the source unattended while the spray voltage is on.

6

Operating in Offline Nanospray Mode

Use offline NSI mode with the Nanospray Flex Series ion source when your LC/MS setup does not include an active sample flow from an LC instrument or the MS syringe pump. For offline NSI mode, mount the offline nanospray source head to the ion source and use Thermo Scientific borosilicate emitters.

Contents

- Offline Nanospray Source Head
- Sample Purification
- Attaching the Offline Nanospray Source Head
- Loading Sample onto Borosilicate Emitters
- Installing the Borosilicate Emitter into the Source Head
- Opening and Positioning the Borosilicate Emitter
- Obtaining a Stable, Low Flow Spray

Offline Nanospray Source Head

The offline nanospray source head (Figure 18), which you can order separately, includes the necessary items for offline analysis. This ion source head contains a custom-made conductive ferrule that makes establishing a proper electrical contact easier between the source head and the borosilicate emitter. A fingertight fitting makes tightening and locking the emitter in the source head also easier. The red PEEK fitting with the 1/16 in. Tefzel[™] tubing provides a fixed gas backpressure.





Sample Purification

The presence of salts, non-volatile buffers, and other polymeric materials (such as those from electrophoresis gels) can affect sample analysis. High amounts of salts might block the emitter, which causes spray instability and low sensitivity due to the presence of salt adducts. This optional purification step can eliminate these issues and concentrate the analytes.

Thermo Fisher Scientific provides a range of StageTips for micro-purification and desalting. For recommended part numbers, see Consumables. The recommended StageTips are based on an Eppendorf[™] GEloader[™] pipette tip with a long narrow tip. To ensure minimal sample loss, you can insert the tip into the borosilicate emitter to directly deposit the sample onto the emitter.

Attaching the Offline Nanospray Source Head



CAUTION Avoid contact with high voltage. Before installing the offline NSI source head, pull the source bottom away from the MS. Then, disconnect the HV cable from the UHPLC liquid junction cross (if installed), and remove the DirectJunction adapter from the XYZ-manipulator arm.

✤ To attach the offline source head to the ion source

Using a 5 mm hex key, remove the hex screw from the **offline nanospray source head**, and then connect the source head to the XYZ-manipulator arm with that screw.



To keep pressure on the syringe, place it in the **static air pressure device** with the backing plate pressed into one of the slots.

Loading Sample onto Borosilicate Emitters

You can load sample onto borosilicate glass emitters either directly or after a purification step. The tools for this procedure are in the Offline Nanospray (Nano ES) Source Head Kit ES259.

* To load sample onto a borosilicate emitter

- 1. Using the precision tweezers, remove an emitter from the transport container.
- 2. Using the glass cutter, shorten the emitter length to minimize contact between the sample and the glass walls.
- 3. (Optional) Use a StageTip to elute the sample directly into the emitter as follows:
 - a. Insert the StageTip pipette tip as far as possible into the emitter without blocking the passage of air that must exit the emitter.

Figure 19. Inserting the StageTip into the borosilicate emitter



b. If the sample is above the tapered end, shake down the liquid by flicking your wrist or use a table centrifuge.

Note Always point the emitter tip downwards to prevent sample from moving up the sides.

Installing the Borosilicate Emitter into the Source Head



CAUTION Avoid personal injury. Before you remove the emitter, make sure that you depressurize the system and disconnect the HV cable to the source bottom. Otherwise, the emitter might eject at a high speed and cause personal injury to you or someone nearby.

To install the borosilicate emitter

Using the precision tweezers, insert the sample-loaded emitter into the

fitting on the source head.





IMPORTANT The internal ferrule is electrically conductive, providing high voltage to the emitter.



Opening and Positioning the Borosilicate Emitter

Thermo Fisher Scientific recommends that you practice the following procedure with standard sample before using more important samples. If you bend the emitter too far, it might break, resulting in an orifice diameter that is too wide and a sample flow that is too high.

* To open and position the borosilicate emitter

- 1. Ensure that the MS is in off mode.
- 2. Open the emitter tip as follows:
 - a. Apply pressure to the sample by pressing the syringe plunger about halfway in.
 - b. Using the XYZ-manipulator arm, offset the emitter from the entrance into the MS.
 - c. Slowly push the source bottom toward the MS until the emitter tip gently touches the MS spray cone (Figure 20).

The slim end of the emitter bends slightly and a small droplet appears when the emitter opens. You can retract this droplet into the emitter by relaxing the air pressure with the syringe. If the emitter does not open, slide the tip vertically or horizontally over the sides of the spray cone.





- 3. Position the emitter tip as follows:
 - a. Using the XYZ-manipulator arm, position the emitter tip 1–3 mm in front of the entrance to the MS spray cone.
 - b. Using your fingers, loosen the black screw on the source head, raise the back of the emitter to an angle of 0–90 degrees above the tip, and then tighten the screw.

Obtaining a Stable, Low Flow Spray

With the syringe in the static air pressure device, apply enough air pressure to obtain a stable signal with a good signal-to-noise ratio. The gas-tight syringe and the spray head create a loss-free system (unless O-rings or ferrules are damaged).

A stable spray typically lasts at least 30 minutes with 1 μ L of liquid. If the spray stops before depleting the sample, the most likely cause is a blocked emitter.

✤ To restart the stopped spray

Try the following:

- a. With the spray voltage turned on, release and reapply the backing air pressure to create a negative backpressure a few times. If there is no effect, briefly apply the maximum backing pressure.
- b. Repeat step a to open the emitter tip but with the spray voltage turned off.

6 Operating in Offline Nanospray Mode Obtaining a Stable, Low Flow Spray

Configuring the Mass Spectrometer for NSI Mode

Follow the applicable procedure in this chapter to configure the Thermo Scientific MS for nanoelectrospray ionization (nanoESI or NSI) mode.

IMPORTANT Before you begin NSI analysis, check that the source's bottom is pushed completely toward the instrument; otherwise, an error message appears.

Contents

- Configuring the NSI Parameters for the Nanospray Flex NG Source
- Configuring the NSI Parameters for the Nanospray Flex Source
- Selecting the Source for the LCQ Deca XP Max Mass Spectrometer

Configuring the NSI Parameters for the Nanospray Flex NG Source

After you complete the instrument configuration, use the Thermo Tune application to configure the NSI source parameters. For additional information, refer to the Tune Help.

- To set the NSI source parameters in Tune
- Open the Tune application by choosing Start > All Apps (Windows 10) or All Programs (Windows 7) > Thermo Instruments (or Thermo *Model*) and so on.
- 2. Click Ion Source, and then click the Ion Source tab to view the Ion Source page.

The MS automatically detects and enters the source type (NSI) in the Ion Source Type box.

3. In the Pos Ion Spray Voltage (V) box, enter 1900.

Use 1900 V as the start value for the spray voltage. If the intensity of the full-scan spectrum is low, gradually increase the spray voltage to improve the spectrum. The recommended range for the spray voltage is 1500–2200 V.

- 4. In the Sweep Gas (Arb) box, enter 0.
- 5. Click Apply.

Configuring the NSI Parameters for the Nanospray Flex Source

After you complete the instrument configuration, use the Thermo Tune Plus application to configure the NSI source parameters. For additional information, refer to the Tune Plus Help.

- ***** To set the NSI source parameters in Tune Plus
- Open the Tune Plus application by choosing Start > All Apps (Windows 10) or All Programs (Windows 7) > Thermo Instruments (or Thermo *Model*) and so on.
- 2. Choose **Setup > NSI Source** to open the NSI Source dialog box.
- 3. In the Spray Voltage (kV) box, enter 1.90.

Use 1.90 kV as the start value for the spray voltage. If the intensity of the full-scan spectrum is low, gradually increase the spray voltage to improve the spectrum. The recommended range for the spray voltage is 1.50–2.20 kV.

4. Click OK.

Selecting the Source for the LCQ Deca XP Max Mass Spectrometer

Note This section is for the LCQ Deca XP Max[™] MS, which uses the Xcalibur[™] data system version 2.0.7 or earlier.

After you install the Nanospray Flex ion source, use the Instrument Configuration window to configure the MS for NSI mode.

To configure the MS for NSI mode

- On the Windows taskbar, choose Start > All Programs > Xcalibur > Instrument Configuration to open the Instrument Configuration window.
- 2. Select the devices to control from the Xcalibur data system if they are not already selected:
 - a. In the Device Types list, select All.
 - b. Under Available Devices, double-click the icons for the MS and nanoelectrospray LC instrument to add them to the Configured Devices list.
- 3. Double-click the instrument icon to open the Model Configuration dialog box.
- 4. In the left pane, select **Ion Source** to display the source configuration page, and then select **Nanospray** in the Default Source list.
- 5. Click OK, and then click OK again to close the message box.
- 6. Configure the LC device if you have not done so.

For instructions, refer to the LC instrument documentation.

- 7. In the Instrument Configuration window, click **Done**.
- 8. Restart the data system computer and the MS.

7 Configuring the Mass Spectrometer for NSI Mode Selecting the Source for the LCQ Deca XP Max Mass Spectrometer

8

Troubleshooting

Typical problems with the Nanospray Flex Series ion source relate to one or both cameras or to the signal intensity. For troubleshooting tips, see Table 5. If you need further assistance, contact your local Thermo Fisher Scientific service engineer. See Contacting Us.

Table 5.	Nanosprav	y Flex	Series id	on source	problems,	causes,	and	possible solutions	(Sheet 1	l of 3	3)
----------	-----------	--------	-----------	-----------	-----------	---------	-----	--------------------	----------	--------	----

Problem Cause		Solution		
Camera issues				
The light is not on. –or– The video output is too dark.	One or both camera lights are off.	Turn on one or both cameras by pressing their light buttons (Figure 3).		
Signal issues				
The signal intensity is weak.	There is a leak somewhere in the liquid path.	Check all of the LC connections.		
	The parameters for the LC, MS, or both might need adjusting.	 Try these solutions: Verify that the emitter is correctly positioned. Check the LC method and MS Tune method parameters. Run a known standard to check the sensitivity. 		
	The MS's spray cone and ion transfer tube are dirty.	Clean both the spray cone and ion transfer tube. For instructions, refer to the instrument documentation.		

Problem	Cause	Solution
The spray is unstable. –or– There is no spray.	There is a leak somewhere in the liquid path.	Check all of the LC connections.
	Air bubbles in the emitter might cause the spray to "spit."	Try degassing the mobile phase or purging the line, and then recheck the line for air bubbles.
	There is an emitter blockage from particles in the sample, other small particles from the flow lines or valves, and so on.	Try adjusting the spray voltage. If that does not resolve the blockage problem, replace the emitter or column.
	The cleaning solvent is not LC/MS grade, which can negatively affect the spray stability.	Use an LC/MS-grade solvent to clean the emitter tip.
	The emitter is out of alignment, which might occur if you bumped the source or moved the source from one MS to another MS.	Follow the procedureTo adjust the emitter tip position.
	The HV connection to the stainless steel emitter might be loose or dirty.	After you place the MS in off mode, check the HV cable connection on the DirectJunction adapter.

 Table 5.
 Nanospray Flex Series ion source problems, causes, and possible solutions (Sheet 2 of 3)

Problem	Cause	Solution	
The spray is unstable. –or– There is no spray.	The source's HV connection might be unstable.	Check the high voltage contact on the back of the sour (Figure 6) and the front of the MS.	
		If the problem continues for the Nanospray Flex NG source, do the following:	
		 Open Tune and view the Ion Source – Ion Source page. 	
		2. Compare the setting for the spray voltage to the adjacent readback value.	
		A green box () indicates that the parameter is functioning properly.	
		If the problem continues for the Nanospray Flex source, do the following:	
		1 2 3	 Open Tune Plus, choose View > Display Status View, and then click the All tab.
			 Choose Setup > NSI Source to open the NSI Source dialog box.
		A green check mark (\checkmark [Spray Voltage (kV)]:) indicates that the parameter is functioning properly.	

 Table 5.
 Nanospray Flex Series ion source problems, causes, and possible solutions (Sheet 3 of 3)

8 Troubleshooting

9

Replaceable Parts

To order any of these parts for the Nanospray Flex Series ion source, contact your local Thermo Fisher Scientific service engineer.

Spare Parts

Camera, analog (composite video) interface, and power supply	ES216
Camera, USB interface, and software disc	ES218
LCD monitor and power supply	ES217
Offline Nano ES ion source head	ES260
Replacement Ferrules and O-Ring Kit	ES294
USB extension cord 00302-99	-00095

Conductive Emitter Configurations

DirectJunction adapter (includes ZDV union SC600)	ES256
Liquid junction Tee 1/32	ES258
Offline Nanospray (Nano ES) Source Head Kit	ES259
Emitters, borosilicate (100 pieces)	ES380
Offline nanospray ion source head	ES260
Precision tweezers and ceramic glass cutter	ES250
Static air pressure device	ES242
Spare parts, including conductive ferrules and O-ring	N/A

Nonconductive Emitter Configurations

HPLC liquid junction cross 1/32	ES257
HPLC micro union kit 1/32, PEEK, red	G10-0035
UHPLC Liquid Junction Kit (360 µm OD tubing)	ES269
Tightening tool	P-278 (IDEX [™])
Tubing, column-out	
Tubing, waste-in	6041.5289
UHPLC blind ferrule plugs	P-116 (IDEX)
UHPLC cross, 360 μm	UH-752 (IDEX)
UHPLC fused-silica union (use with ES269)	ES272
UHPLC microferrules	PK-152 (IDEX)

Consumables

Emitters, SilicaTip [™] , uncoated FS360-20-10-N-5-105CT (New Ob	jective)
Emitters, borosilicate, long (1.1 mm tip), 52 mm long, Au/Pd double coating,	
(100 pieces)	ES388
Emitters, borosilicate, medium (0.8 mm tip), 52 mm long, Au/Pd double coating,	
(100 pieces)	ES387
Emitters, borosilicate, medium (0.8 mm tip), 52 mm long, Au/Pd single coating,	
(100 pieces)	ES380
StageTips, C8, GELoader pipette tip (96 pieces)	SP121
StageTips, C18, GELoader pipette tip (96 pieces)	SP101
StageTips, SCX, GELoader pipette tip (96 pieces)	SP141
For other consumables, visit www.fishersci.com and www.unitylabservices.com.	

A

Contents of the Installation Kits

Table 6 lists the parts supplied in the Nanospray Flex NG Ion Source Kit (P/N ES072) and the Nanospray Flex Ion Source Kit (P/N ES071), which are identical except for the source housing. For a list of replaceable parts, see Chapter 9, "Replaceable Parts."

Note As of 2018, these ion source kits exclude the LCD monitor (P/N ES217) and the 1.5 mm hex key.

Image	Part	Quantity	Part number
Provided in the Nanospray Flex NG Ion Source Kit (P/N ES072)			
See Figure 2.	Nanospray Flex NG housing	1	_a
Provided in the Nanospray	Flex Ion Source Kit (P/N ES071)		
See Figure 3.	Nanospray Flex housing	1	_b
Provided in both ion source	e kits		
	Dino-Lite analog cameras and power supplies	2 ^c	ES216 or
	–or– Dino-Lite USB cameras and software disc; USB extension cords		ES218 (cameras); 00302-99-00095 (cords)
	DirectJunction Adapter	1	ES256
	UHPLC Liquid Junction Kit ^d	1	ES269

Table 6. Contents of the Nanospray Flex Series ion source kits (Sheet 1 of 2)

Image	Part	Quantity	Part number
_	Liquid Junction Column Out Line (tubing), 360 µm OD, 20 µm ID, 550 mm long	1	6041.5290
-	Liquid Junction Waste In Line (tubing), 360 µm OD, 75 µm ID, 550 mm long	1	6041.5289
	Emitter, stainless steel, 150 µm OD, 30 µm ID, 40 mm long (4 pieces)	1	ES542
_	Emitters, New Objective SilicaTip, uncoated, 360 µm OD, 20 µm ID, 10.5 cm long (5 pieces)	1	FS360-20- 10N5- 105CT
-	Emitter sleeves, PEEK, 1/32 in. OD, 280 µm (10 pieces)	1	SC903
-	Hex key, 3 mm	1	-
-	Spare parts (blind plug, O-rings, screws, sleeves)	_	-

Table 6. Contents of the Nanospray Flex Series ion source kits (Sheet 2 of 2)

^a The Nanospray Flex NG source housing comes with the Nanospray Flex NG Ion Source Kit (P/N ES072) and cannot be purchased separately.

^b The Nanospray Flex source housing comes with the Nanospray Flex Ion Source Kit (P/N ES071) and cannot be purchased separately.

^c The ion source kit includes either two of the analog cameras (P/N ES216) with one LCD monitor (P/N ES217) or two of the USB cameras with two USB extension cords.

^d In 2015, the UHPLC kit replaced the Liquid Junction Cross 1/32 (P/N ES257).

Index

A

analysis, offline 2

C

cameras DinoCapture video application (USB camera) 10 focusing wheel 11 mounting bracket (USB camera) 10 troubleshooting 41 type provided 9 *See also* procedures column, configuring one or two 14 compliance regulatory xiii consumables, part numbers 45 contacting us xiv

D

DinoCapture application (USB camera) 12 directive, WEEE iii DirectJunction adapter column configuration 14 *See also* procedures documentation online survey x documentation, accessing ix downloading documents x

E

emitters borosilicate (offline mode) 32 setup angle and distance 26 stainless steel or glass 13 *See also* procedures

F

figures, list of vii

G

gas connection, optional 8

H

high voltage caution statement 28 connector, source back 7 DirectJunction cable 27 troubleshooting 42

ion sources compatible mass spectrometers xi locking levers, positions 6 optional gas connection 8 photo of 3–4, 7 *See also* procedures

L

Liquid Junction Cross 1/32 48

Μ

mass spectrometer ion sweep cone, types 6 NSI mode, configuring 37 preparing before installing the source 5 spray cone attached to MS 7

Ν

nanoelectrospray, advantages 2 NSI, offline flow rate 2 NSI, online flow rate 2

0

offline analysis 2 offline mode 29

P

procedures cameras, installing 9 DirectJunction adapter, attaching 25 emitters borosilicate installing 33 opening and positioning the tip 34 glass (HPLC), configuring 19 glass (UHPLC), configuring 20 glass or stainless steel tip position, adjusting 27 offline borosilicate, loading sample 32 stainless steel (HPLC), configuring 16-17 ion source attaching to the MS 6 removing from the MS 8 monitor, installing 9 nanoViper fitting, using 15 nitrogen gas (optional), connecting 8 offline source head, installing 31 offline spray, restarting 35 video picture, adjusting 12

R

regulatory compliance xiii replaceable parts 45 resources, additional 23

S

safety high voltage 28 symbol descriptions xi signal issues, troubleshooting 41 source *See* ion sources spray voltage, setting the value 38 symbols, meaning xi

Т

Thermo Scientific website, user documents x troubleshooting 10, 41 Tune application, opening 37 Tune Plus (legacy) application, opening 38

U

UHPLC liquid junction cross, mounting 22

W

WEEE directive iii

X

XYZ manipulator control knobs 28