

H-ESI Probe

User Guide

97055-97045 Revision C April 2009

© 2009 Thermo Fisher Scientific Inc. All rights reserved.

The following are registered trademarks in the United States and other countries:

Teflon and Tefzel are registered trademarks of E. I. du Pont de Nemours & Co. KEL-F is a registered trademark of 3M Corporation.

PEEK is a trademark of Victrex plc.

Ion Max is a trademark of Thermo Fisher Scientific Inc. in the United States. Xcalibur is a registered trademark of Thermo Fisher Scientific Inc. in the United States.

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, June 2005; Revision B, March 2008; Revision C, April 2009

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

WEEE Compliance

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



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

WEEE Konformität

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



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

Contents

	Preface	vii
	Safety and Special Notices	vii
	Precautions for Handling the H-ESI Probe	viii
	Contacting Us	ix
Chapter 1	Introduction	1
Chapter 2	Functional Description	5
Chapter 3	Removing and Installing the H-ESI Probe	7
	Removing the H-ESI Probe	7
	Installing the H-ESI Probe	9
Chapter 4	Maintenance	13
	Flushing the Sample Transfer Line, Sample Tube, and H-ESI Probe	13
	Trimming the H-ESI Sample Tube	14
	Replacing the H-ESI Needle and Needle Seal	14
	Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve	16
Chapter 5	Replaceable Parts	19
	Index	21

Preface

This *H-ESI Probe User Guide* provides you with information on using the heated-electrospray ionization (H-ESI) probe. It also provides procedures for installing and maintaining the H-ESI probe.



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

Safety and Special Notices

Make sure you follow the precautionary statements presented in this guide. The safety and other special notices appear in boxes.

Safety and special notices include the following:



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

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

Note Highlights information of general interest.

Tip Highlights helpful information that can make a task easier.

Precautions for Handling the H-ESI Probe

When operating the mass spectrometer in the heated-electrospray mode, do **not** touch the heated surfaces of the H-ESI probe or the Ion Max API source housing. Touching the heated surfaces of the probe or the source housing can cause burns when the probe is operating at typical temperatures of 350 to 450 °C.



CAUTION AVOID BURNS. At typical operating temperatures (350 to 450 °C), the H-ESI probe and the API source housing can severely burn you. Before removing the probe from the API source housing, allow the probe to cool to room temperature (for approximately 20 minutes) before you touch it.

When operating the mass spectrometer in the heated-electrospray mode, if you observe liquid leaking from the sample inlet, place the mass spectrometer in Standby mode before you tighten the fitting to eliminate the leak.



CAUTION AVOID ELECTRIC SHOCK. Do not tighten the probe sample inlet fitting to eliminate a liquid leak while the mass spectrometer is in operation. If you touch liquid leaking from the probe sample inlet while the mass spectrometer is in operation, you might receive an electric shock.

Contacting Us

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

❖ To contact Technical Support

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

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

❖ To contact Customer Service for ordering information

Phone	800-532-4752
Fax	561-688-8731
E-mail	us.customer-support.analyze@thermofisher.com
Web site	www.thermo.com/ms

❖ To copy manuals from the Internet

Go to mssupport.thermo.com and click **Customer Manuals** in the left margin of the window.

❖ To suggest changes to documentation or to Help

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



- Send an e-mail message to the Technical Publications Editor at techpubs-lcms@thermofisher.com.

Introduction

Heated-electrospray ionization (H-ESI) transforms ions in solution into ions in the gas phase by using electrospray ionization (ESI) in combination with heated auxiliary gas. H-ESI can be used to analyze any polar compound that makes a preformed ion in solution. The term *preformed ion* can include adduct ions. For example, polyethylene glycols can be analyzed from a solution containing ammonium acetate because of adduct formation between the NH_4^+ ions in the solution and oxygen atoms in the polymer.

This chapter describes the principles of H-ESI and the H-ESI probe. See [Figure 1](#). The H-ESI probe is housed in the Ion Max™ ion source housing.

Figure 1. Ion source interface (left) and H-ESI probe (right)



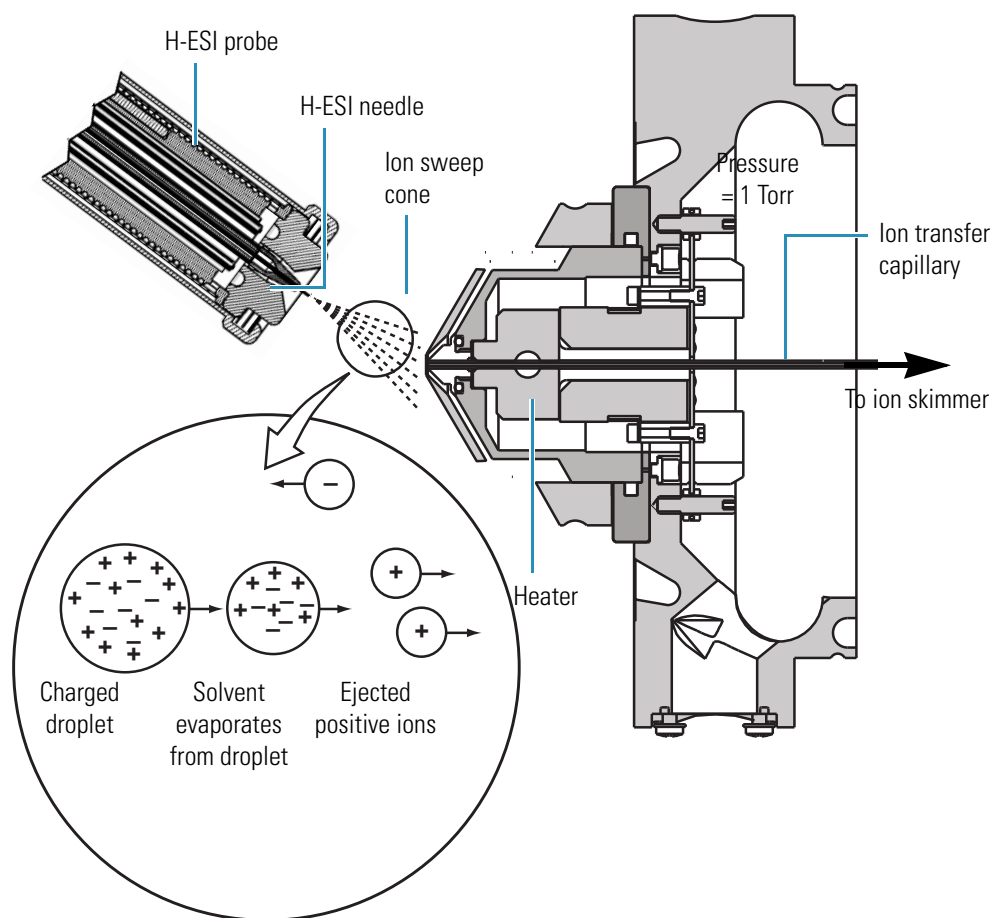
With H-ESI, a mass spectrometer can analyze a range of molecular weights greater than 100 000 u, due to multiple charging. H-ESI is especially useful for the mass analysis of polar compounds, which include biological polymers (for example, proteins, peptides, glycoproteins, and nucleotides); pharmaceuticals and their metabolites; and industrial polymers (for example, polyethylene glycols).

In H-ESI, ions are produced and analyzed as follows:

1. The sample solution enters the H-ESI needle, which receives a high voltage.
2. The H-ESI needle sprays the sample solution into a fine mist of droplets that are electrically charged at their surface.
3. The electrical charge density at the surface of the droplets increases as solvent evaporates from the droplets. In H-ESI, heated auxiliary gas aids solvent evaporation.
4. The electrical charge density at the surface of the droplets increases to a critical point known as the Rayleigh stability limit. At this critical point, the droplets divide into smaller droplets because the electrostatic repulsion is greater than the surface tension. The process is repeated many times to form very small droplets.
5. Electrostatic repulsion ejects sample ions from the very small, highly charged droplets into the gas phase.
6. The sample ions enter the mass spectrometer and are analyzed.

Figure 2 shows the steps in the formation of ions from highly charged droplets.

Figure 2. H-ESI process in the positive ion polarity mode



You can use the H-ESI mode in either positive or negative ion polarity mode. The ion polarity mode of choice is determined by the polarity of the preformed ions in solution: Acidic molecules form negative ions in solution, and basic molecules form positive ions. Use a positively charged needle to analyze positive ions and a negatively charged needle to analyze negative ions, as the ejection of sample ions from droplets is facilitated if the ionic charge and surface charge of the droplet have the same polarity.

Sample ions can carry a single charge or multiple charges. The number of charges carried by the sample ion depends on the structure of the analyte of interest and the carrier solvent. (In H-ESI, the buffer and the buffer strength both have a noticeable effect on sensitivity, so it is important to choose these variables correctly.) In the case of higher molecular weight proteins or peptides, the resulting mass spectrum consists typically of a series of peaks corresponding to a distribution of multiply charged analyte ions.

Droplet size, surface charge, liquid surface tension, solvent volatility, and ion solvation strength are factors that affect the H-ESI process. Large droplets with high surface tension, low volatility, strong ion solvation, low surface charge, and high conductivity prevent good electrospray.

Organic solvents, such as methanol, acetonitrile, and isopropyl alcohol, are superior to water for H-ESI. Volatile acids and bases are good, but salts above 10 mM concentration and strong acids and bases are extremely detrimental.

Follow these rules for achieving a good electrospray:

- Keep salts out of the solvent system.
- Use organic/aqueous solvent systems and volatile acids and bases.
- Optimize the pH of the solvent system.

Note You do not need to tune and calibrate the mass spectrometer as part of your daily routine.

Calibration parameters are instrument parameters that affect the mass accuracy and resolution. Tune parameters are instrument parameters that affect the intensity of the ion signal. You must tune and calibrate the mass spectrometer (that is, optimize the tune and calibration parameters) approximately once a quarter. Refer to the “Automatic Tuning and Calibrating in the H-ESI/MS Mode” chapter in your instrument’s getting started manual for a procedure to tune and calibrate your mass spectrometer.

You must optimize the tune parameters (or change the Tune Method) whenever you change the type of experiment. Refer to the “Optimizing the Mass Spectrometer with Your Compound in H-ESI/MS Mode” chapter in your instrument’s getting started manual for a procedure to optimize and tune the parameters for your H-ESI experiment.

Table 1 shows initial H-ESI settings for different liquid flow rates. These initial settings provide a starting point for optimizing system performance. The optimal settings for your application depend on the compounds of interest, the solvent matrix, and the chromatographic conditions. For information on optimizing these settings, refer to the getting started guide for your mass spectrometer.

For an LTQ Series mass spectrometer, the allowable range for the auxiliary gas flow depends on the version of the instrument control software and the vaporizer temperature setting:

- For LTQ 2.5.0 or lower, the allowable range for the auxiliary gas flow is 5 to 60 units.
- For LTQ 2.5.5 or higher, the allowable range for the auxiliary gas flow is 0 to 60 units for vaporizer temperatures up to 100.00 °C. For vaporizer temperatures above 100.00 °C, the minimum auxiliary gas flow is 5 units.

For a TSQ Series mass spectrometer, you can set the auxiliary gas flow from 0 to 60 units.

Tip For best results, set the auxiliary gas flow to a minimum of 5 units for vaporizer temperatures above 100.00 °C.

Table 1. Initial H-ESI settings for the H-ESI probe

Liquid flow rate (µL/min)	Ion transfer tube (capillary) temperature (°C)*	H-ESI vaporizer temperature (°C)**	Sheath gas pressure (psi)	Auxiliary gas flow (arbitrary units)	Spray voltage (V)
5	240	Off to 50	5	0 ^{***}	+3000 (-2500) ^{****}
200	350	250 to 300	35	30	+3000 (-2500)
500	380	300 to 400	60	50	+3000 (-2500)
1000	400	350 to 450	75	60	+3000 (-2500)

* Always optimize the tube lens voltage whenever you change the temperature of the ion transfer tube.

** Compound dependent

*** For an LTQ Series mass spectrometer controlled from LTQ 2.5.0 or lower, the minimum auxiliary gas flow setting is 5.

**** Negative ion mode

Functional Description

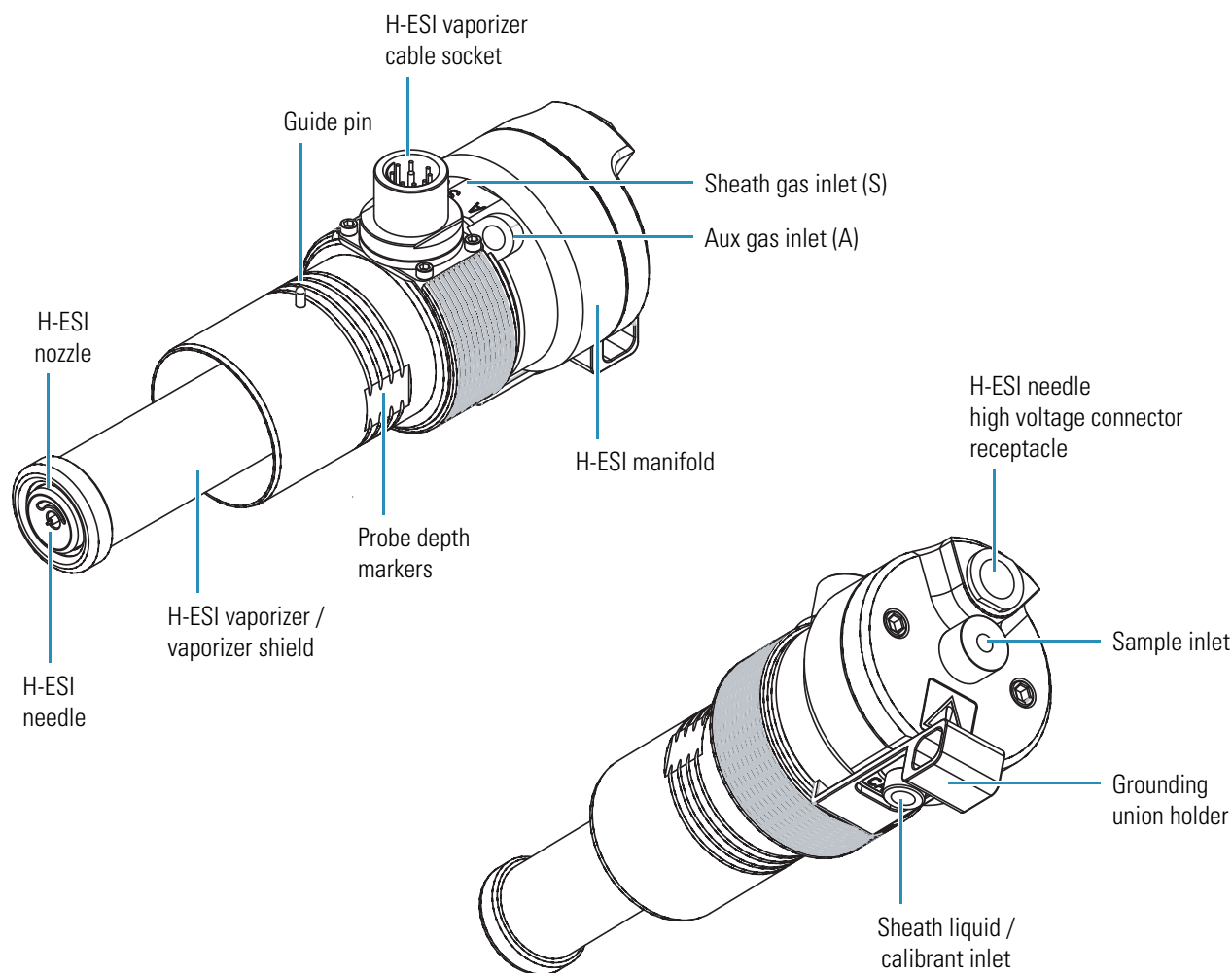
The H-ESI probe, shown in [Figure 3](#), produces charged aerosol droplets that contain sample ions. The H-ESI probe accommodates liquid flows of 1 $\mu\text{L}/\text{min}$ to 1 mL/min without splitting.

The H-ESI probe includes a fused-silica sample tube, needle, nozzle, manifold, vaporizer, and heat exchanger. Sample and solvent enter the H-ESI probe through the sample tube. The sample tube is a short section of 0.1-mm ID, fused-silica tubing that extends through the H-ESI probe and into the H-ESI needle to near the end of the H-ESI needle. A large negative or positive voltage is applied to the H-ESI needle (typically ± 3 to ± 5 kV), which sprays the sample solution into a fine mist of charged droplets. The H-ESI nozzle directs the flow of sheath gas and auxiliary gas at the droplets. The H-ESI manifold houses the H-ESI nozzle and needle, and includes the sheath gas and auxiliary gas plumbing. The sheath gas plumbing delivers dry nitrogen gas to spray the sample solution out the nozzle.

The H-ESI probe has inlets for the introduction of sample solution, sheath gas, auxiliary gas, and sheath liquid or calibrant into the H-ESI probe. The sheath gas is the inner coaxial nitrogen gas that sprays (nebulizes) the sample solution into a fine mist as it exits the sample tube. Typical sheath gas flow rates for H-ESI are 5 to 20 units for sample flow rates of less than 10 $\mu\text{L}/\text{min}$, and 50 to 75 units for sample flow rates greater than 400 $\mu\text{L}/\text{min}$. When you tune the mass spectrometer, for best results adjust the sheath gas flow rate until the ion signal is stable.

The heated auxiliary gas is the outer coaxial nitrogen gas that assists the sheath gas in the desolvation of sample solutions. The auxiliary gas heats as it spirals through a vaporizer. The vaporizer is thermally insulated from the sample tube to prevent direct heating of the sample solution. You can control the temperature of the vaporizer in the Xcalibur™ data system. The temperature range is from room temperature to 600 °C. For recommended operating temperatures and gas flow settings, see [Table 1](#) on [page 4](#).

Figure 3. H-ESI probe



To enhance desolvation or to improve the stability of the electrospray, you can infuse sheath liquid directly from the syringe pump into the H-ESI probe via the sheath liquid/calibrant inlet (labeled CAL). You can also infuse internal calibrants into the H-ESI probe.

The angle of the H-ESI probe is fixed at approximately 60 degrees. To help optimize the spray stability, use adjustment screws on the Ion Max source housing to make small changes to probe position. The fixed angle, off-axis spraying affords long-term signal stability (robustness) for most solutions that contain non-volatile matrix components, mobile phase buffers, or ion-pairing reagents. For information on adjusting the probe position, refer to the *Ion Max and Ion Max-S API Source Hardware Manual*.

Removing and Installing the H-ESI Probe

This chapter describes how to remove and install the H-ESI probe from the Ion Max source housing. You must remove the H-ESI probe to change ionization modes and to perform maintenance on the probe. You do not need to remove the Ion Max source housing to change probes.

Note To remove or install an APCI or ESI probe, refer to the *Ion Max and Ion Max-S API Source Hardware Manual*.

Removing the H-ESI Probe



CAUTION AVOID BURNS. At operating temperatures, the H-ESI vaporizer can severely burn you. The H-ESI vaporizer typically operates between 350 and 450 °C. **Always allow the heated vaporizer to cool to room temperature (for approximately 20 min) before you remove or touch the H-ESI probe.**

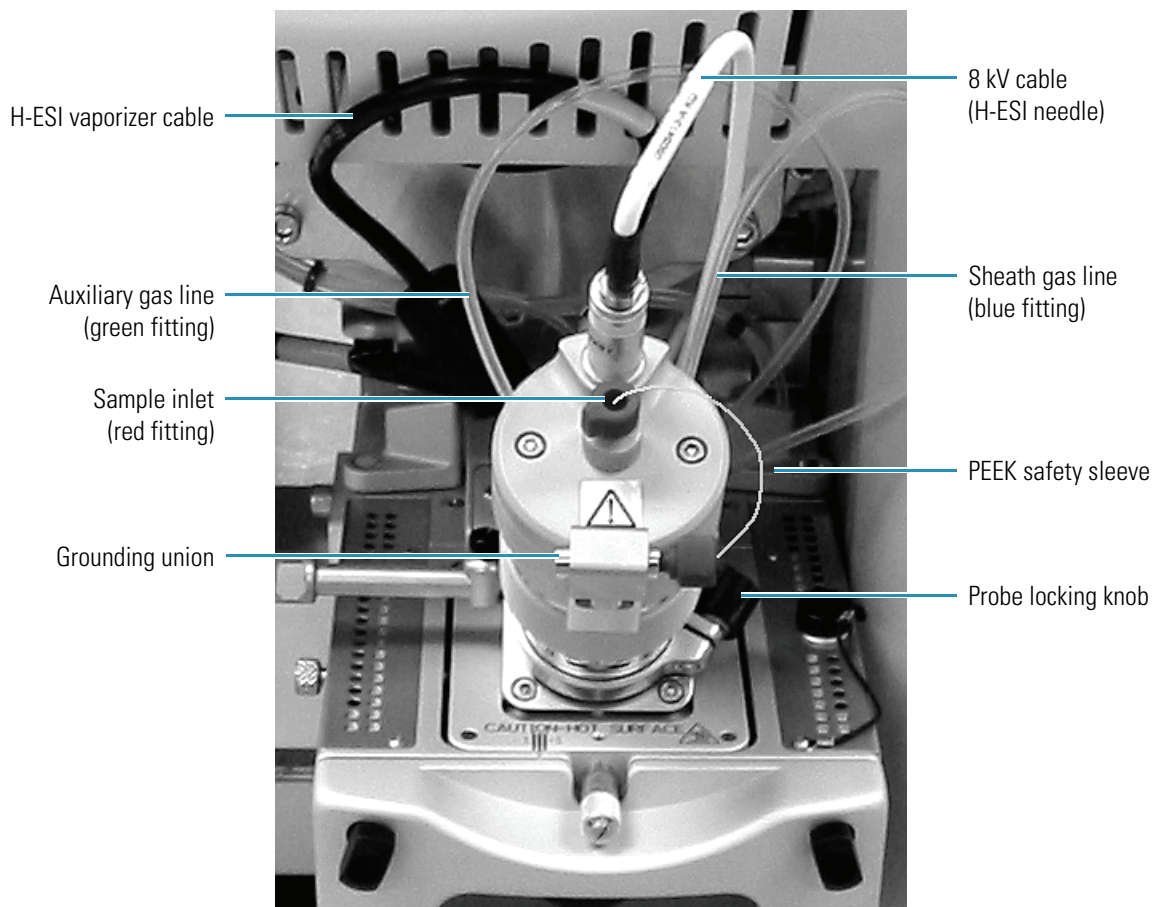
❖ To remove the H-ESI probe

1. Place your mass spectrometer in standby.
2. Disconnect the sample transfer tubing (LC line) from the grounding union (stainless steel ZDV fitting).
3. Disconnect the 8 kV cable from the H-ESI needle high voltage receptacle as follows (see [Figure 4](#)):
 - a. Unlock the cable by twisting the locking ring counterclockwise.
 - b. Unplug the 8 kV cable from the H-ESI needle high voltage receptacle.
4. Unplug the H-ESI vaporizer cable from the cable socket on the H-ESI probe. See [Figure 4](#).
5. Connect the H-ESI vaporizer cable to the ESI interlock socket on the ion source housing. See [Figure 5](#).
6. Disconnect the auxiliary gas fitting (green) from the auxiliary gas inlet (A) on the probe manifold. See [Figure 4](#).
7. Disconnect the sheath gas fitting (blue) from the sheath gas inlet (S) on the probe manifold.

3 Removing and Installing the H-ESI Probe

Removing the H-ESI Probe

Figure 4. H-ESI probe installed on the Ion Max housing



8. Unlock the probe locking ring by turning the probe locking knob counterclockwise.
9. Carefully pull the probe straight back in the port in the housing until it meets with the slot in the ESI interlock block. The guide pin on the probe manifold prevents you from twisting the probe until the pin is aligned with the slot in the ESI interlock block. Once the probe is all the way back and aligned with the slot, turn the probe 45 degrees counterclockwise to free the probe from the alignment notch. Be careful not to break the fused-silica sample tube or PEEK™ safety sleeve.
10. Pull the probe straight out to remove it from the ion source housing.
11. Store the H-ESI probe in its original shipping container.

Installing the H-ESI Probe

Note To remove an APCI or ESI probe, refer to the *Ion Max and Ion Max-S API Source Hardware Manual*.

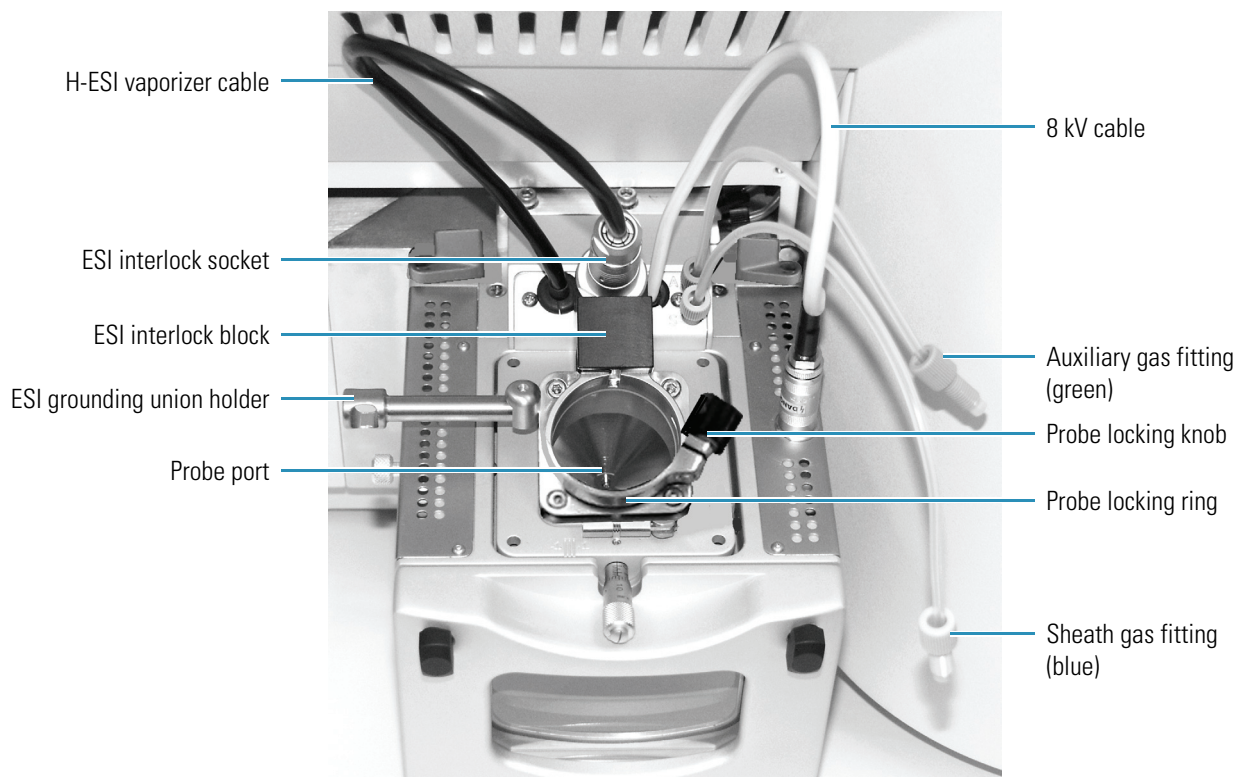
❖ **To install the H-ESI probe**

1. Remove the H-ESI probe from its storage container. Inspect and clean it if necessary.

Note If your H-ESI probe does not already have a sample tube (fused-silica capillary) and safety sleeve attached, you must follow the procedure for installing a sample tube and PEEK safety sleeve in the topic “[Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve](#)” on page 16.

2. Ensure that the probe locking ring is opened to its widest position. See [Figure 5](#).

Figure 5. Ion Max ion source housing without H-ESI probe

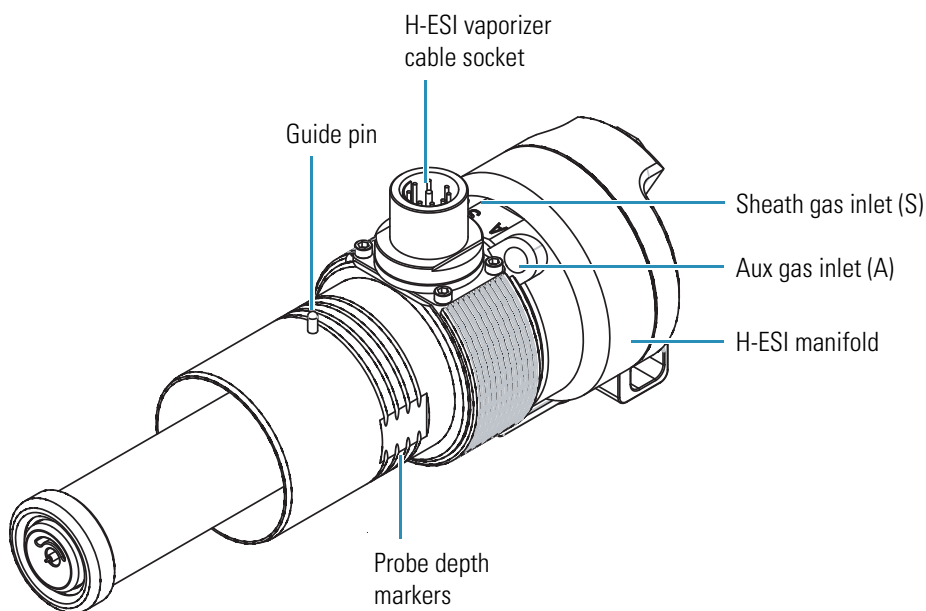


3 Removing and Installing the H-ESI Probe

Installing the H-ESI Probe

3. Insert the H-ESI probe into the port in the ion source housing. Make sure to align the guide pin on the probe body (see [Figure 6](#)) at a minus 45° angle from the ESI interlock block (see [Figure 5](#)).

Figure 6. H-ESI probe, front view

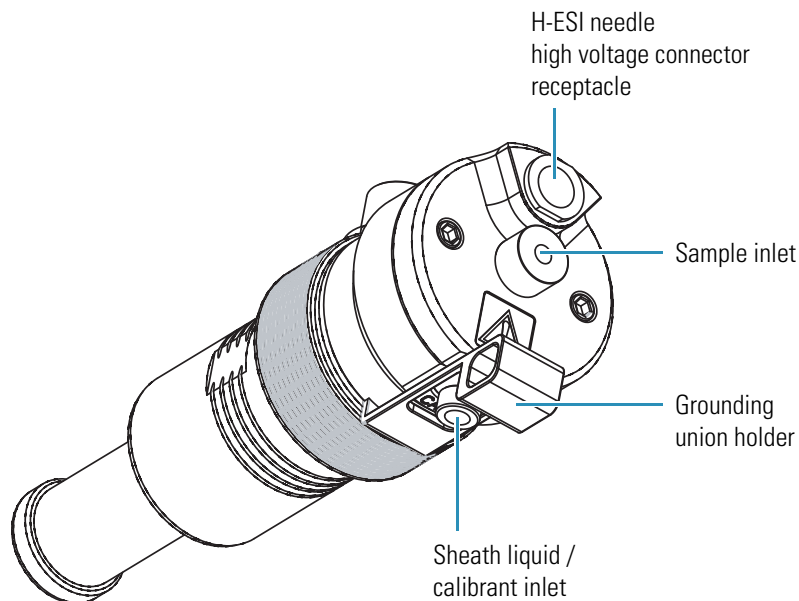


4. Push the probe into the port until the guide pin meets with the locking ring on the ion source housing. See [Figure 5](#).
5. Turn the probe 45 degrees clockwise and align the guide pin with the slot in the ESI interlock block; you might have to pull the probe towards you slightly to properly align the pin with the notch. Once you have turned the probe far enough to align the pin with the alignment notch at the rear of the port, push the probe straight in until the guide pin stops at the bottom of the alignment notch.
6. Lock the probe in place by turning the probe locking knob clockwise. See [Figure 5](#).
7. Ensure that the grounding union (stainless steel ZDV fitting) is seated in the grounding union holder on the H-ESI probe. See [Figure 7](#).



CAUTION To avoid electric shock, ensure that the grounding union is seated in the grounding union holder on the H-ESI probe. If the grounding union is not seated and you touch the fused-silica solvent line while the system is in operation, you might receive an electric shock.

Figure 7. H-ESI probe, rear view

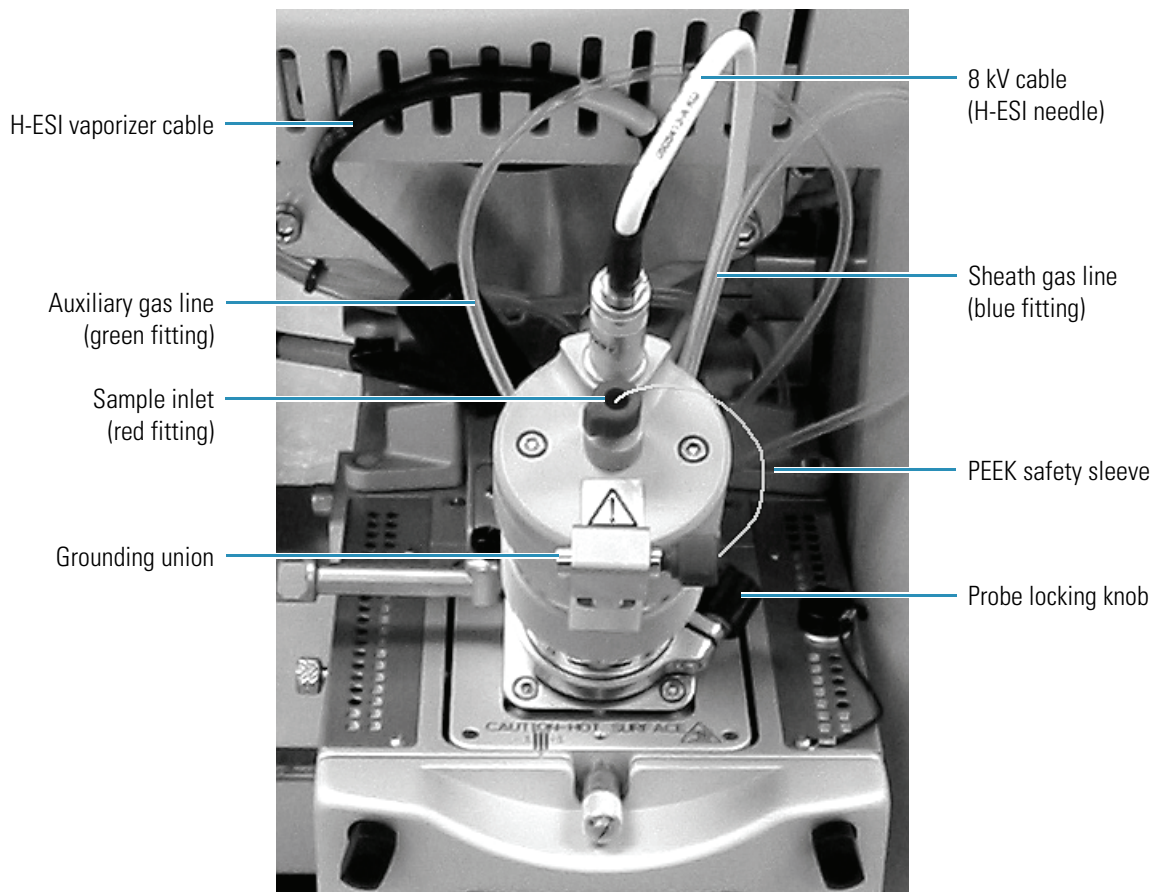


8. Connect the sheath gas fitting (blue) to the sheath gas inlet (S) on the probe manifold. See [Figure 6](#).
9. Connect the auxiliary gas fitting (green) to the auxiliary gas inlet (A) on the probe manifold. See [Figure 6](#).
10. Unplug the H-ESI vaporizer cable from the ESI interlock socket.
11. Connect the H-ESI vaporizer cable to the vaporizer cable socket on the H-ESI probe. See [Figure 6](#).
12. Connect the 8 kV cable to the H-ESI needle high voltage receptacle on the H-ESI probe. See [Figure 7](#). Tighten the locking ring on the 8 kV connector.
13. Connect the sample transfer tubing (LC line) to the grounding union. See [Figure 8](#).

3 Removing and Installing the H-ESI Probe

Installing the H-ESI Probe

Figure 8. H-ESI probe installed in the Ion Max housing



The H-ESI source is now properly installed on the mass spectrometer.

Tip Before you analyze samples with the H-ESI source, you must change to H-ESI source mode in Quantum Tune Master by choosing **Setup > Change Ion Source > HESI**.

Maintenance

The H-ESI probe requires minimum maintenance. If the fused-silica sample tube is plugged or broken, you must replace it. Although you can trim or replace the sample tube without disassembling the H-ESI probe, to replace the H-ESI needle or needle seal, you must partially disassemble the H-ESI probe.

Tip For best results, flush the H-ESI probe at the end of each working day, using a 50:50 LCMS-grade methanol\distilled water solution from the LC through the H-ESI probe.

Wear clean gloves when you handle H-ESI probe components.



CAUTION AVOID BURNS. At operating temperatures, the H-ESI vaporizer can severely burn you. The H-ESI vaporizer typically operates between 350 and 450 °C. Always allow the heated vaporizer to cool to room temperature (for approximately 20 min) before you remove or touch the H-ESI probe.

Contents

- [Flushing the Sample Transfer Line, Sample Tube, and H-ESI Probe](#)
- [Trimming the H-ESI Sample Tube](#)
- [Replacing the H-ESI Needle and Needle Seal](#)
- [Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve](#)

Flushing the Sample Transfer Line, Sample Tube, and H-ESI Probe

For best results flush the sample transfer line, sample tube, and H-ESI probe for 15 minutes at the end of each working day (or more often if you suspect they are contaminated). Use a 50:50 methanol\distilled water solution from the LC through the API source. After 15 minutes, turn off the flow of liquid from the LC to the API source, but keep the API source on (including the sheath gas and auxiliary gas) for an additional 5 minutes. Refer to the daily operations chapter in your mass spectrometer's hardware manual.

Trimming the H-ESI Sample Tube

Operating your instrument with acetonitrile in the mobile phase can cause the polyimide coating on the fused-silica sample tube to elongate. If the polyimide coating has elongated past the end of the H-ESI needle, you must cut and reposition the end of the sample tube.

❖ **To cut and reposition the end of the sample tube approximately 0.5 mm (between 0 and 1 mm) inside the end of the H-ESI needle**

1. Remove the H-ESI probe from the Ion Max source by following the procedure in the topic [“Removing the H-ESI Probe”](#) on [page 7](#).
2. Loosen the sample inlet fitting.
3. Gently pull back on the sample tube to free it from the fitting.
4. Push the sample tube forward so that it extends beyond the end of the H-ESI needle.
5. Use a fused-silica cutting tool to cut off a small length of sample tube. Ensure that you cut the end of the sample tube squarely.
6. Pull the sample tube backwards until the exit end of the sample tube is recessed just inside the H-ESI needle by approximately 0.5 mm (between 0 and 1 mm).
7. Tighten the sample inlet fitting securely to hold the sample tube.

Note The sample tube might move forward when you tighten the sample inlet fitting. Ensure that the sample tube is retracted into the H-ESI needle approximately 0.5 mm (between 0 and 1 mm). If necessary, loosen the fitting and reposition the sample tube.

8. Reinstall the H-ESI probe as described in the topic [“Installing the H-ESI Probe”](#) on [page 9](#).

Replacing the H-ESI Needle and Needle Seal

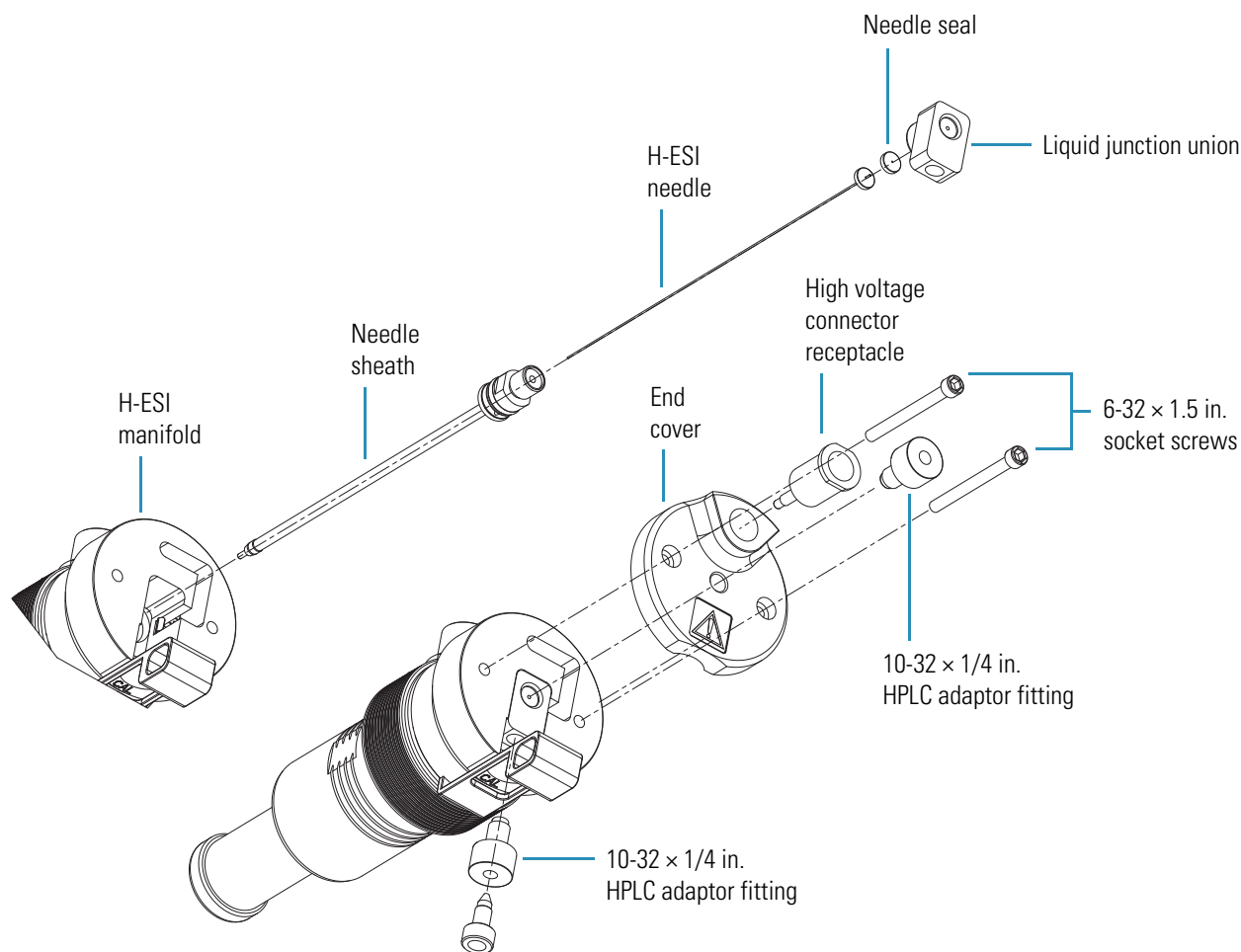
You must replace the H-ESI needle if it is damaged. You need to replace the needle seal if the sheath gas is leaking at the needle seal-needle interface. If you replace the needle, you should also replace the needle seal.

❖ **To replace the needle and/or needle seal (see [Figure 9](#))**

1. Remove the H-ESI probe from the Ion Max source by following the procedure in the topic [“Removing the H-ESI Probe”](#) on [page 7](#).
2. Unscrew the sample inlet fitting.
3. Remove the sample tube and sample inlet fitting from the H-ESI probe.
4. Use a 7/64 in. hex wrench or ball driver to remove the two 6-32 × 1/4 in. socket screws.
5. Remove the end cover from the H-ESI manifold.

6. Unscrew and remove the sheath liquid/calibrant fitting.
7. Remove the liquid junction union, needle, needle seal, and needle sheath from the H-ESI manifold as a unit. See [Figure 9](#).
8. While grasping the liquid junction union with a 1/2 in. open end wrench, use a 3/8 in. open end wrench to unscrew the needle sheath from the liquid junction union.
9. Remove the H-ESI needle and needle seal from the needle sheath.
10. Remove the needle seal from the H-ESI needle.
11. Inspect the Teflon™ needle seal (P/N 97055-20271). If the needle seal is deformed, replace it.
12. Inspect the 26-gauge H-ESI needle (P/N 97055-20273). If the H-ESI needle is damaged, replace it. If you replace the needle, you should also replace the needle seal.
13. Reinstall the H-ESI probe as described in the topic [“Installing the H-ESI Probe”](#) on [page 9](#).

Figure 9. H-ESI probe (exploded view)



❖ To reassemble the H-ESI probe (see [Figure 9](#))

1. Insert the entrance end of the H-ESI needle into the needle seal.
2. Reassemble the needle sheath, H-ESI needle and seal, and liquid union junction unit as follows:
 - a. Seat the H-ESI needle and needle seal in the needle sheath.
 - b. Thread the needle sheath into the liquid junction union. Slightly wet the needle sheath threads with HPLC-grade methanol for lubrication.
 - c. With a 3/8 in. wrench, gently tighten the needle sheath until it is a little more than fingertight. Do not overtighten the needle sheath.
3. Insert the needle sheath, H-ESI needle and seal, and liquid union junction unit into the H-ESI manifold.
4. Insert the sheath liquid/calibrant adaptor into the H-ESI manifold and tighten until fingertight.
5. Position the end cover on the H-ESI manifold.
6. Insert the two 6-32 × 1/4 in. socket screws into the H-ESI probe and tighten them with a 7/64 in. hex wrench or ball driver.
7. Carefully insert the fused-silica sample tube into the H-ESI probe.
8. Tighten the sample inlet fitting.
9. Reinstall the H-ESI probe as described in the topic [“Installing the H-ESI Probe”](#) on [page 9](#).

Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve

❖ To install a new sample tube and PEEK safety sleeve

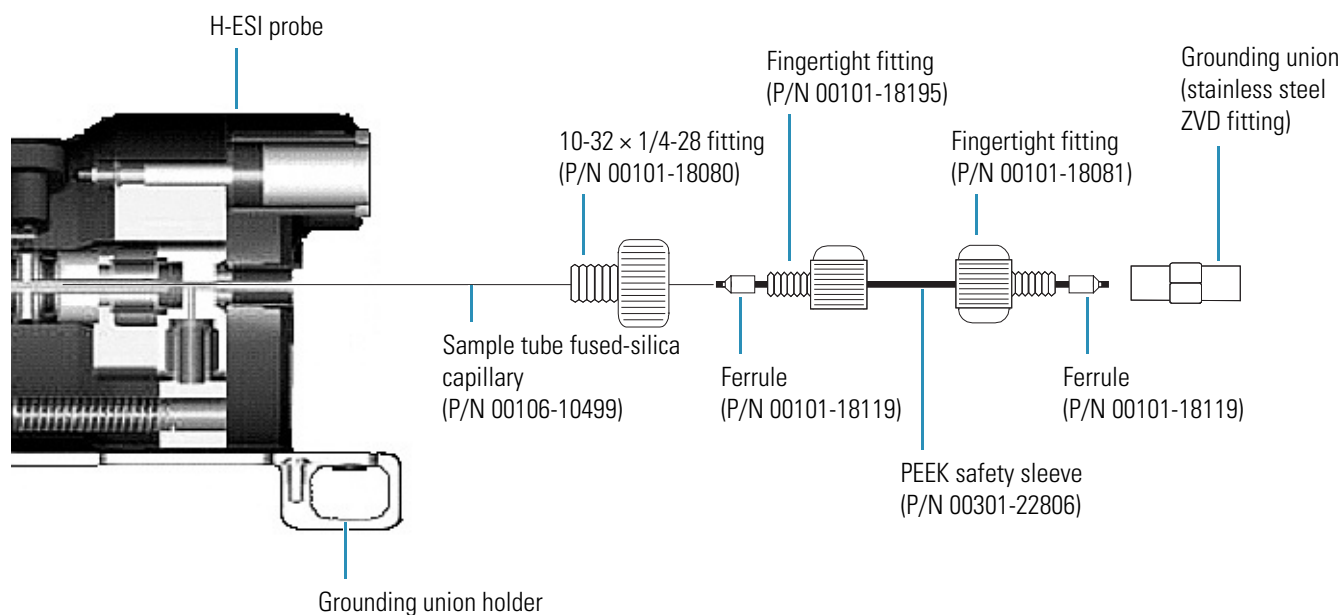


CAUTION AVOID ELECTRICAL SHOCK. When you are operating your instrument in the H-ESI mode, you could receive an electrical shock unless you install the safety kit discussed below. You could receive an electrical shock if the fused-silica capillary tube breaks during H-ESI operation. Therefore, for your safety and in compliance with international safety standards, you **must** cover the fused-silica capillary tube with the PEEK safety sleeve (P/N 00301-22806) and associated PEEK ferrules (P/N 00101-18119) provided in the Safety Sleeve Kit (P/N 70005-62015) before you operate the instrument. Installation instructions (P/N 70005-97009) are included in the kit. Operation of the instrument without the safety sleeve impairs the safety protection provided by the instrument and, thus, could lead to serious injury.

1. Remove the H-ESI probe from the Ion Max source by following the procedure in the topic [“Removing the H-ESI Probe”](#) on [page 7](#).

- Use a fused-silica cutting tool to cut a 30 cm (12 in.) piece of 0.1 mm ID \times 0.19 mm OD, fused-silica tubing (sample tube) (P/N 00106-10499). Ensure that you cut the ends of the fused-silica tubing squarely.
- Insert the sample tube through the exit end of the H-ESI needle and into the H-ESI probe.
- Push the sample tube through the H-ESI probe until approximately 3.5 cm (1.5 in.) is left protruding from the exit end of the H-ESI needle. The remaining length of sample tube should exit the H-ESI probe sample inlet.
- Slide the 10-32 \times 1/4-28 Kel-F™ fitting adaptor (P/N 00101-18080) over the sample tube (see Figure 10) and tighten the fitting onto the H-ESI probe sample inlet.

Figure 10. Fused-silica sample tube and safety sleeve assembly



- Slide the precut 25 cm (10 in.) 0.009 in. ID \times 0.024 in. OD PEEK safety sleeve (P/N 00301-22806) over the sample tube.
- Slide the 0.027 in. ID PEEK ferrule (P/N 00101-18119), narrow end first, over the PEEK safety sleeve and to the 10-32 \times 1/4-28 Kel-F fitting adaptor.
- Slide the (red) Fingertight fitting (P/N 00101-18195) onto the PEEK safety sleeve and into the H-ESI probe sample inlet (labeled *Sample*). Tighten the fitting slightly, but not completely.
- Push the PEEK safety sleeve over the sample tube until it stops against the Teflon needle seal inside the H-ESI probe.
- Pull the sample tube (from the H-ESI needle end) until the sample tube is flush with the precut square end of the PEEK safety sleeve.

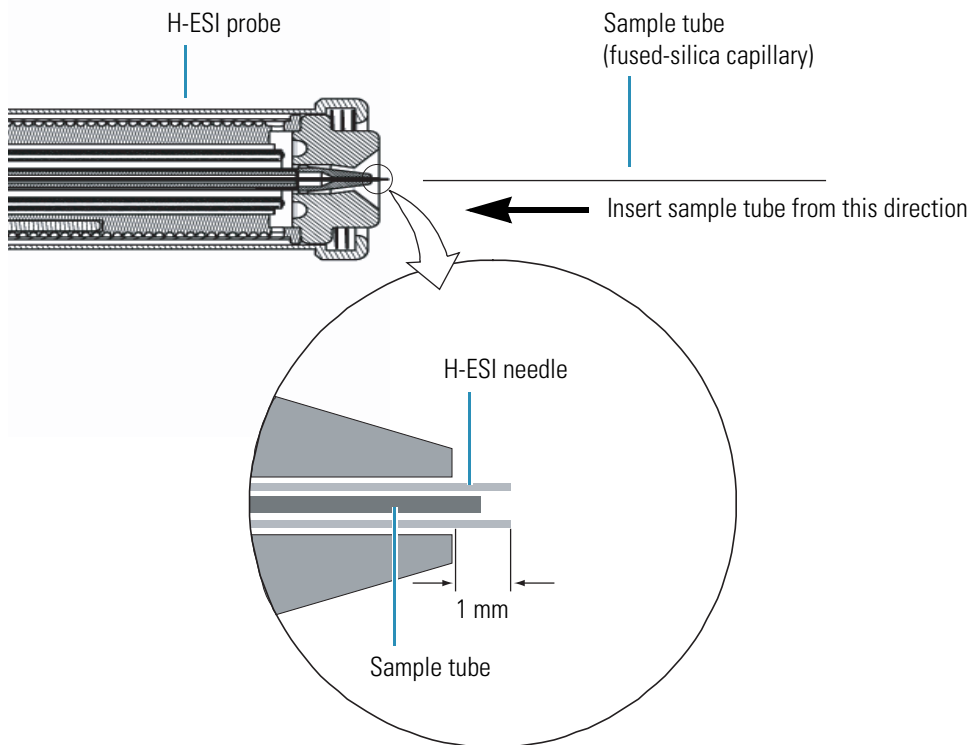
4 Maintenance

Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve

11. Slide a (brown) Fingertight fitting (P/N 00101-18081) and (brown) ferrule (P/N 00101-18119), wide end first, over the PEEK safety sleeve.
12. Connect the PEEK safety sleeve and the ferrule to the (stainless steel) ZDV fitting by tightening the (brown) Fingertight fitting. Ensure that the Fingertight fitting is securely tightened around the PEEK safety sleeve; otherwise, the sample stream might enter between the sample tube and the PEEK safety sleeve. Ensure the sample tube is held tightly in the grounded fitting by gently pulling the sample tube from the exit end of the H-ESI needle.
13. Use a fused-silica cutting tool to cut the sample tube at the H-ESI needle so that only 2.5 cm (1 in.) remains protruding from the exit end of the H-ESI needle.
14. From the sample inlet pull the PEEK safety sleeve backwards, so that the exit end of the sample tube is recessed just inside the H-ESI needle by approximately 0.5 mm (between 0 and 1 mm). See [Figure 11](#).
15. Tighten the (red) Fingertight fitting securely to hold the PEEK safety sleeve and sample tube in place.

Note The sample tube might move forward when you tighten the sample inlet fitting. Ensure that the sample tube is retracted into the H-ESI needle approximately 0.5 mm (between 0 and 1 mm). If necessary, loosen the fitting and reposition the sample tube.

Figure 11. Installing the fused-silica sample tube



Replaceable Parts

This chapter contains part numbers for replaceable and consumable parts for the H-ESI probe. To ensure proper results in servicing the H-ESI probe, order only the parts listed or their equivalent.

This chapter contains part numbers for the following items:

Probe, H-ESI	97055-60140
Needle, metal, 26-gauge, H-ESI	97055-20273
Seal, needle	97055-20271
Tubing, fused-silica, 0.10 mm ID × 0.19 mm OD	00106-10499
Fitting, grounding union, ZDV, 1/4-28	00109-00304
Ferrule, 0.008 in. ID, KEL-F, HPLC	00101-18075
Fitting, plug, 1/4-28, TEFZEL™, HPLC	00101-18075
Ferrule, 0.012 in.ID, KEL-F, HPLC	00101-18116
Safety Sleeve Kit	70005-62015
Ferrule 0.027 in. ID PEEK HPLC	00101-18119
Safety sleeve 0.009 in. ID × 0.024 in. OD × 10 in. length, natural PEEK	00301-22806
Instructions safety sleeve	70005-97009
Fitting, Fingertight Upchurch	00101-18195
Fitting, Fingertight, HPLC, 10-32, PEEK	00101-18081
Fitting, adaptor, 10-32 × 1/4-28, KEL-F	00101-18080

For information on ordering LCMS-grade solvents from Thermo Fisher Scientific, go to www.FisherLCMS.com.

Figure 12. Consumables for the H-ESI probe

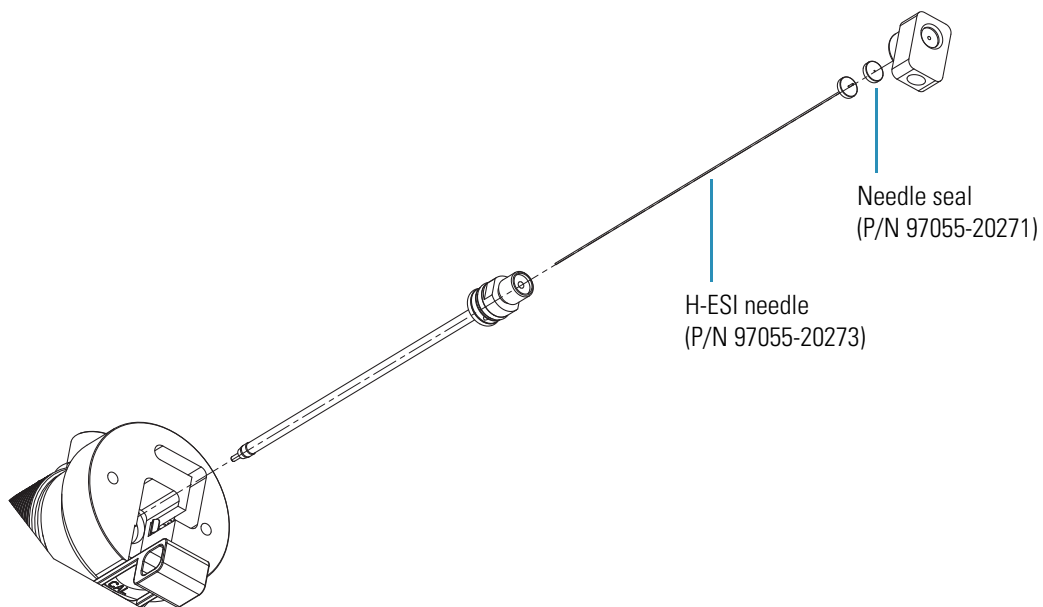
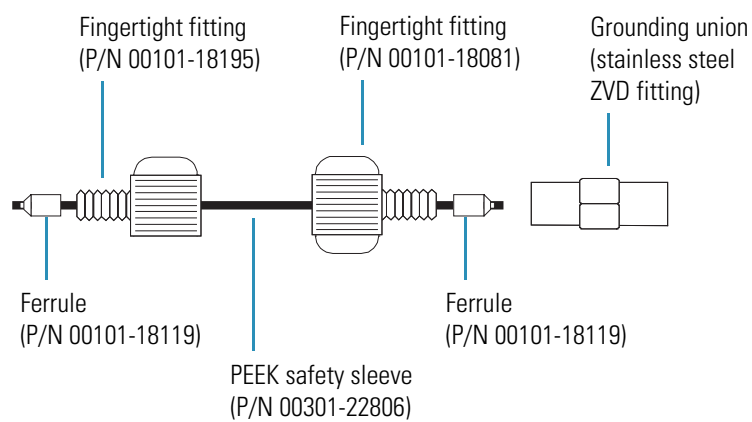


Figure 13. PEEK safety sleeve part numbers



Index

A

APCI probe, removing (note) 7
 auxiliary gas, flow rate settings 4, 5

C

calibration frequency (note) 3
 CAUTIONS
 avoiding an electric shock viii
 avoiding burns, H-ESI vaporizer 13
 electrical shock, PEEK safety sleeve 16
 cleaning procedures
 H-ESI probe, flushing 13
 sample transfer line, flushing 13
 sample tube, flushing 13
 compliance
 WEEE iii

D

daily maintenance
 H-ESI probe, flushing 13
 sample transfer line, flushing 13
 sample tube, flushing 13

E

electrospray, rules for good electrospray 3
 ESI probe, removing (note) 7

F

figures
 H-ESI probe 1, 6
 H-ESI probe exploded view 15
 H-ESI process 2
 installing PEEK safety sleeve 17
 installing sample tube 18
 ion source interface 1
 fused-silica sample tube
 installing 16
 installing (figure) 17, 18
 trimming 14

G

guide pin location (figure) 6

H

heated auxiliary gas, description 5
 H-ESI
 change ion source (tip) 12
 process (figure) 2
 process, discussed 1
 rules for optimizing 3
 H-ESI needle
 high voltage connector, location (figure) 6
 replacing 14
 voltages 5
 H-ESI probe
 disassembling 14
 exploded view (figure) 15
 figure 1, 10
 flushing 13
 functional description 5
 installed (figure) 8, 12
 installing 9
 maintenance overview 13
 PEEK safety sleeve 16
 reassembling 16
 removing 7
 replaceable parts 19
 sample tube, description of 5
 sample tube, installing 16
 H-ESI sample tube, trimming 14
 H-ESI vaporizer
 avoiding burns (CAUTION) 13
 cable socket, location (figure) 6

I

Ion Max (figure) 8, 9, 12
 ion source interface (figure) 1

M

maintenance

- H-ESI probe, overview 13
- needle seal, replacing 14
- needle, replacing 14
- sample transfer line, flushing 13
- sample tube, flushing 13
- sample tube, installing 16
- sample tube, trimming 14

N

needle seal

- location (figure) 15
- reinstalling 16
- removing 14

needle, H-ESI

- location (figure) 15
- reinstalling 16
- removing 14

P

part numbers for replaceable parts 19

PEEK safety sleeve

- CAUTION 16
- installing 16
- installing (figure) 17

R

replaceable parts for the H-ESI probe 19

S

safety precautions for handling the HESI-II probe viii

sample inlet, location of (figure) 6

sample transfer line, flushing 13

sample tube

- flushing 13
- installing 16
- installing (figure) 18
- repositioning (note) 18
- trimming 14

sheath gas

- description 5
- flow rates 5

sheath liquid, description of 6

solvents, ordering 19

T

tuning frequency (note) 3

V

voltages applied to the H-ESI needle 5

W

WEEE compliance iii