LTQ XL ®

Getting Started

97355-97042 Revision A

June 2006



For Research Use Only Not for use in Diagnostic Procedures



© 2004 Thermo Electron Corporation. All rights reserved.

Microsoft[°] and Windows[°] are registered trademarks of Microsoft Corporation. Teflon[°] is a registered trademark of E.I. du Pont de Nemours & Company. Unimetrics[°] is a registered trademark of Unimetrics Company. Tygon[°] is a registered trade mark of Saint-Gobain Performance Plastics Company. Rheodyne[°] and the Rheodyne logo (see) are registered trademarks of Rheodyne, L.P.

All other trademarks are the property of Thermo Electron Corporation and its subsidiaries.

Technical information contained in this publication is for reference purposes only and is subject to change without notice. Every effort has been made to supply complete and accurate information; however, Thermo Electron Corporation assumes no responsibility and will not be liable for any errors, omissions, damage, or loss that might result from any use of this manual or the information contained therein (even if this information is properly followed and problems still arise).

This publication is not part of the Agreement of Sale between Thermo Electron Corporation and the purchaser of an LC/MS system. In the event of any conflict between the provisions of this document and those contained in Thermo Electron Corporation's Terms and Conditions, the provisions of the Terms and Conditions shall govern.

System Configurations and Specifications supersede all previous information and are subject to change without notice.

Printing History: Revision A printed June 2006. Software Revision: LTQ 2.2, Xcalibur 2.2



Regulatory Compliance

Thermo Electron San Jose performs complete testing and evaluation of its products to ensure full compliance with applicable domestic and international regulations. When the system is delivered to you, it meets all pertinent electromagnetic compatibility (EMC) and safety standards as described below.

EMC Directive 89/336/EEC

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

EN 55011	1999	EN 61000-4-3	2002	EN 55011	1998
		IEC 61000-4-3	A1-1998		
EN 61000-3-2	1995, A1; 1998, A2; 1998, A14; 2000	EN 61000-4-4 IEC 61000-4-4	1995, A1; 2001, A2; 2001; A2-1995	EN 61000-3-2	1995, A1; 1998, A2; 1998, A14; 2000
EN 61000-3-3	1998	EN 61000-4-5	1995, A1; 2001	EN 61000-3-3	1998
		IEC 61000-4-5	2005		
EN 61326-1	1998, A3	EN 61000-4-6	1996, A1; 2001	EN 61326-1	1998
		IEC 61000-4-6	2004		
EN 61000-4-2	2000	EN 61000-4-11	1994, A1; 2001	EN 61000-4-2	2000
IEC 61000-4-2	2001	IEC 61000-4-11	2001-03		
FCC Class A, CFR 47 Part 18	2005	CISPR 11	1999, A1; 1999, A2; 2002		

Low Voltage Safety Compliance

This device complies with Low Voltage Directive EN 61010-1:2001.

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



FCC Compliance Statement

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



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

Notice on Lifting and Handling of Thermo Electron San Jose Instruments

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

Notice on the Proper Use of Thermo Electron San Jose Instruments

In compliance with international regulations: If this instrument is used in a manner not specified by Thermo Electron San Jose, the protection provided by the instrument could be impaired.



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

WEEE Konformität

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



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



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 Electron 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 Electron à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Electron qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur <u>www.thermo.com/WEEERoHS</u>.

Contents

Preface	xi
About This Guide	xi
Related Documentation	xi
Safety and Special Notices	xi
Contacting Us	xii
Assistance	xii
Changes to the Manual and Online Help	xii

Chapter 1	Introduction1
	Why Use the LTQ XL MS Detector?
	Which MS Detector Technique—ESI or APCI—Is Better for
	Analyzing My Samples?4
	Using ESI/MS
	Using APCI/MS
	Should I Use Sheath, Auxiliary, and/or Sweep Gases?7
	How Can I Introduce My Samples into the MS Detector?
	What Types of Buffers Should I Use? What Types Should I
	Avoid?
	How Should I Set Up the MS Detector for Various LC Flow
	Rates?
	What is Tuning and Calibration of the MS Detector All About?
	13
	What Types of Experiments Can I Perform with the LTQ XL MS
	Detector?
	General MS or MSn Experiments16
	Data-Dependent Experiments
	Ion Mapping Experiments
	Ion Tree Experiments
	1

Chapter 2 Setting Up the Ion Source for Tuning and Calibrating the MS Detector27

Placing the LC/MS System in Standby	
Removing the APCI Probe	29
Removing the Ion Max Ion Source Housing	
Installing the Ion Sweep Cone	
Installing the Ion Max Ion Source Housing	
Installing the ESI Probe	
motaning the Let Trobe	

Chapter 3	Tuning and Calibrating Automatically in the ESI/MS Mode 43 Setting Up the Syringe Pump for Tuning and Calibration
	Cleaning the MS Detector after Tuning and Calibrating62
Chapter 4	Tuning with Your Analyte in LC/ESI/MS Mode69Setting Up to Introduce Sample by Syringe Pump into Solvent Flow from an LC71Setting Up to Tune the MS Detector with Your Analyte74Optimizing the MS Detector Tune Automatically with Your Analyte77Saving the ESI/MS Tune Method81
Chapter 5	Acquiring ESI Sample Data Using the Tune Plus Window83 Setting Up to Acquire MS/MS Data in the Full Scan Type84 Optimizing the Isolation Width and Setting Up to Optimize the Collision Energy
Chapter 6	Setting Up the Ion Source for Acquiring Data in APCI/MS/MS Mode101Removing the ESI Probe102Removing the Ion Max Ion Source Housing105Removing the Ion Sweep Cone106Installing the Corona Needle107Installing the Ion Max Ion Source Housing108Installing the APCI Probe111
Chapter 7	Optimizing the MS Detector with Your Analyte in APCI/MS Mode115 Setting Up the Inlet for Tuning Using High-Flow Infusion116 Setting Up the MS Detector for APCI/MS Operation

	Optimizing the Tune of the MS Detector Automatically in APCI/MS Mode
Chapter 8	Acquiring APCI Sample Data Using the Tune Plus Window131
	Setting Up to Introduce Sample by Loop Injection into Solvent
	Flow from an LC132
	Acquiring APCI Data in the SIM Scan Mode134
Appendix A	Sample Formulations139
	Caffeine, MRFA, and Ultramark 1621 Stock Solutions
	Stock Solution: Caffeine
	Stock Solution: MRFA142
	Stock Solution: Ultramark 1621142
	ESI Calibration Solution: Caffeine, MRFA, Ultramark 1621143
	Reserpine
	Stock Solution: Reserpine144
	Reserpine Tuning Solution and Reserpine APCI Sample
	Solution144
Appendix B	LTQ XL High Mass Range Calibration147
	High Mass Range Calibration Solution
	Normal Mass Range Calibration
	Enter Normal Mass Range Data into Tune Plus
	Tune on m/z 524.3
	High Mass Range Calibration Procedure154
	Notes

Preface

About This Guide	Welcome to the Thermo Electron, LTQ XL™ system. The LTQ XL is a member of the Thermo family of MS detectors.		
	This <i>LTQ XL Getting Started</i> manual provides information on how to set up calibrate, and tune the LTQ XL MS detector, and how to acquire LC/MS data. All of these procedures can be performed from the Xcalibur [®] Tune Plu window.		
	To perform analyses in ESI mode, see Chapters 2, 3, 4, and 5. To perform analyses in APCI mode, see Chapters 2, 3, 6, 7, and 8.		
Related Documentation	In addition to this guide, Thermo Electron provides the following documents for the LTQ XL system:		
	LTQ XL Preinstallation Guide		
	LTQ XL Getting Connected		
	• LTQ XL Hardware Manual		
	Help is also available from within the software.		
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 laser-related hazards to human beings. It includes information specific to the class of laser involved. Each DANGER notice is accompanied by the international laser radiation symbol.		
	CAUTION Highlights begands to humans, property, or the opving mont		
	Each CAUTION notice is accompanied by an appropriate CAUTION symbol.		

	IMPORTANT Highlights information necessary to avoid damage to software, loss of data, invalid test results, or information critical for optimal performance of the system.			
	Note Highlights information of general interest.			
	Note Helpful information that can make a task easier.			
Contacting Us	There are several ways to contact Thermo Electron.			
Assistance	For new product updates, technical support, and ordering information, contact us in one of the following ways:			
	Visit Us on the Web			
	www.thermo.com/finnigan			
	Contact Technical Support			
	Phone: 1-800-685-9535 Fax: 1-561-688-8736 techsupport.finnigan@thermo.com			
	Contact Customer Service			
	In the US and Canada for ordering information:Phone:1-800-532-4752Fax:1-561-688-8731			
	International contacts for ordering information: Visit www.thermo.com/finnigan for the current listing,			
Changes to the Manual and Online Help	To suggest changes to this guide or to the Help, use either of the following methods:			
	• Fill out a reader survey online at www.thermo.com/lcms-techpubs			
	 Send an e-mail message to the Technical Publications Editor at techpubs.finnigan-lcms@thermo.com 			

Chapter 1 Introduction

The LTQ XL[™] is a member of the Thermo family of MS detectors. The LTQ XL MS detector is an advanced analytical instrument that includes a syringe pump, a divert/inject valve, an atmospheric pressure ionization (API) source, an MS detector, and the Xcalibur data system. In a typical analysis, a sample can be introduced in any one of the following ways:

- Using the syringe pump (direct infusion)
- Using the inject valve fitted with a loop and an LC pump (flow injection analysis)
- Using a valve and an LC system fitted with a column (LC/MS)

In analysis by LC/MS, a sample is injected onto an LC column. The sample is then separated into its various components. The components elute from the LC column and pass into the MS detector where they are analyzed. Analysis by direct infusion or flow injection provides no chromatographic separation of components in the sample before it passes into the MS detector. The data from the MS detector is then stored and processed by the Xcalibur data system.

This introduction answers the following questions:

- Why Use the LTQ XL MS Detector?
- Which MS Detector Technique—ESI or APCI—Is Better for Analyzing My Samples?
- How Can I Introduce My Samples into the MS Detector?
- What Types of Buffers Should I Use? What Types Should I Avoid?
- How Should I Set Up the MS Detector for Various LC Flow Rates?
- What is Tuning and Calibration of the MS Detector All About?
- What Types of Experiments Can I Perform with the LTQ XL MS Detector?

Why Use the LTQ XL MS Detector?

The attribute that sets the LTQ XL MS detector apart from other LC detectors is the high level of analytical specificity that it provides. The LTQ XL MS detector can provide multiple levels of analysis. Each level of analysis adds a new dimension of specificity for positive compound identification. The various levels of analysis are as follows:

- Chromatographic separation and compound detection (non MS technique utilizing chromatographic retention time)
- Mass analysis (molecular mass information)
- Two-stage mass analysis, MS/MS (structural information)
- Multiple MS-MS mass analysis, MSⁿ (structural information)
- ZoomScan[™] analysis (charge state information)

Chromatographic separation and compound detection can be obtained by all LC/detector systems. Retention time alone, however, does not positively identify a compound because many compounds can have the same retention time under the same experimental conditions. In addition, even if a compound is identified correctly by retention time, quantitation results can be in error because other compounds in the sample might co-elute with the compound of interest.

Single stage mass analysis allows for the identification of analytes of interest. Atmospheric pressure ionization typically produces mass spectra that provide molecular mass information.

Two-stage mass analysis allows for even more positive compound identification. MS/MS analysis monitors how a parent ion fragments when exposed to an additional stage of ionization. There are two types of MS/MS analysis: Full Scan MS/MS and Selective Reaction Monitoring (SRM). Full Scan MS/MS monitors the production of all product ion from a specific parent ion. SRM MS/MS analysis monitors a specific reaction path: the production of a specific product ion from a specific parent ion. Using MS/MS analysis, you can easily quantitate target analytes in complex matrices such as plant or animal tissue, plasma, urine, groundwater, or soil. Because of the specificity of MS/MS measurements and the ability to eliminate interferences by an initial mass selection stage, quantitative target compound analysis is easily accomplished using the LTQ XL MS detector.

Multiple MS/MS mass analysis provides a unique capability to obtain structural information that can be useful in structure elucidation of metabolites, natural products, and sugars. MSⁿ techniques on the LTQ XL MS detector allow for stepwise fragmentation pathways, making interpretation of MSⁿ spectra relatively straightforward. The LTQ XL MS detector has several advanced features that make its MSⁿ capabilities extremely powerful for qualitative analysis. (See "What Types of Experiments Can I Perform with the LTQ XL MS Detector?" on page 16.)

ZoomScan analysis provides information about the charge state of one or more mass ions of interest. ZoomScan data is collected by using slower scans at higher resolution. This allows for unambiguous determination of charge state, which in turn allows for the correct determination of molecular mass.

In addition to the aforementioned levels of analysis, there is an additional technique called Wideband Activation. The Wideband Activation option allows the LTQ XL MS detector to apply collision energy to ions during MS/MS fragmentation over a fixed mass range of 20 u. This option allows the LTQ XL MS detector to apply collision energy to both the parent ion, as well as to product ions created as a result of non-specific losses of water (18 u) or ammonia (17 u), for example, or to product ions formed from the loss of fragments less than 20 u. When you want enhanced structural information and you do not want to perform MS³ analysis with the LTQ XL MS detector, choose the Wideband Activation option for qualitative MS/MS. Because the collision energy is applied to a broad mass range, signal sensitivity is somewhat reduced when you choose this option. Therefore, increase the value of the collision energy (Activation Amplitude) to compensate somewhat for the reduction of sensitivity.

Which MS Detector Technique—ESI or APCI—Is Better for Analyzing My Samples?

The LTQ XL MS detector includes two standard atmospheric pressure ionization source probes: $^{\rm 1}$

- Electrospray ionization (ESI) probe
- Atmospheric pressure chemical ionization (APCI) probe

Typically, more polar compounds such as amines, peptides, and proteins are best analyzed by ESI, and nonpolar compounds such as steroids are best analyzed by APCI.

Sample ions can carry a single charge or multiple charges. The number of charges carried by the sample ions depends on the structure of the analyte of interest, the mobile phase, and the ionization mode.

Using ESI/MS The ESI mode typically produces mass spectra consisting of multiply charged ions (for proteins and peptides) depending on the structure of the analyte and the solvent. For example, the resulting mass spectrum of a higher molecular mass protein or peptide typically consists of a distribution of multiply charged analyte ions. The resulting mass spectrum can be mathematically manipulated to determine the molecular mass of the sample.

The ESI mode transfers ions in solution into the gas phase. Many samples that previously were not suitable for mass analysis (for example, heat-labile compounds or high molecular mass compounds) can be analyzed by ESI. 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 NH₄⁺ ions in the solution and oxygen atoms in the polymer. With ESI, the range of molecular masses that can be analyzed by the LTQ XL MS detector is greater than 100,000 u, due to multiple charging. 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.

You can use the ESI mode in either positive or negative ion polarity mode. The ion polarity mode is determined by the polarity of the preformed ions in solution: Acidic molecules form negative ions in high pH solution, and basic molecules form positive ions in low pH solution. A positively charged ESI needle is used to generate positive ions and a negatively charged needle is used to generate negative ions.

¹Optional ionization sources [atmospheric photo ionization (APPI), atmospheric pressure matrix assisted laser desorption ionization (AP MALDI), and nanospray] are also available.

You can vary the flow rate from the LC into the MS detector over a range from 1 μ L/min to 1000 μ L/min. See Table 3. (In ESI, the buffer and the buffer strength both have a noticeable effect on sensitivity. Therefore, it is important to choose these variables correctly.) In the case of higher molecular mass proteins or peptides, the resulting mass spectrum consists typically of a series of peaks corresponding to a distribution of multiply charged analyte ions.

The ESI process is affected by droplet size, surface charge, liquid surface tension, solvent volatility, and ion solvation strength. Large droplets with high surface tension, low volatility, strong ion solvation, low surface charge, and high conductivity prevent good electrospray.

Mixed organic/aqueous solvent systems that include organic solvents such as methanol, acetonitrile, and isopropyl alcohol are superior to water alone for ESI. Volatile acids and bases are good, but salts above 10 mM are not recommended. Strong mineral acids and bases are extremely detrimental to the instrument.

The rules for a good electrospray are as follows:

- Keep non-volatile salts and buffers out of the solvent system. For example, avoid the use of salts containing sodium or potassium and avoid the use of phosphates. If necessary, use ammonium salts instead.
- Use organic/aqueous solvent systems and volatile acids and bases.
- If possible, optimize the pH of the solvent system for your analyte of interest. For example, if your analyte of interest contains a primary or secondary amine, your mobile phase should be slightly acidic (pH 2 to 5). The acid pH tends to keep positive ions in solution.

Using APCI/MS Like ESI, APCI is a soft ionization technique. APCI provides molecular mass information for compounds of medium polarity that have some volatility. APCI is typically used to analyze small molecules with molecular masses up to about 2000 Da.

APCI is a gas phase ionization technique. Therefore, the gas phase acidities and basicities of the analyte and solvent vapor play an important role in the APCI process.

APCI is a very robust ionization technique. It is not affected by minor changes in most variables such as changes in buffer or buffer strength. The rate of solvent flowing from the LC into the MS detector in APCI mode is typically high (between 0.2 and 2 mL/min). See Table 3.

You can use APCI in positive or negative ion polarity mode. For most molecules, the positive-ion mode produces a stronger ion current. This is especially true for molecules with one or more basic nitrogen (or other basic) atoms. Molecules which generally produce strong negative ions, with acidic sites such as carboxylic acids and acid alcohols, are an exception to this general rule.

Although, in general, fewer negative ions are produced than positive ions, negative ion polarity can be more specific. This is because the negative ion polarity mode sometimes generates less chemical noise than does the positive mode. Thus, the signal-to-noise ratio might be better in the negative ion mode than in the positive ion mode.

Should I Use Sheath, Auxiliary, and/or Sweep Gases?

Nitrogen gas can be applied to the system using any combination of the three gas sources: Auxiliary gas, Sweep gas, and/or Sheath gas. When Sheath gas is used, nitrogen is applied as an inner coaxial gas (when used in tandem with Auxiliary gas), helping to nebulize the sample solution into a fine mist as the sample solution exits the ESI or APCI nozzle. (Sheath gas is not used with the NSI source.) When Auxiliary gas is being used, nitrogen flows through the ion source nozzle, the vapor plume is affected; the spray is focused and desolvation is improved. When Sweep gas is used, the nitrogen flows out from behind the sweep cone and can result in solvent declustering and adduct reduction.

When you are analyzing complex matrices such as plasma or nonvolatile salt buffers, Sweep gas is required for ruggedness. In full-scan MS or data dependent scan experiments, the signal-to-noise ratio can be improved by application of Sweep gas. In some cases, signal intensity can be increased by using Auxiliary gas, particularly for higher LC flow rates.

All analyses are analyte dependent and require separate optimization with Sheath gas, Sweep gas, and Auxiliary gas to determine which combination will yield optimum performance. It is especially important to optimize with each gas independently before you perform experiments using MSⁿ techniques and before you perform any quantitative analysis experiments because optimum results could be achieved with any combination of Sheath, Sweep, and/or Auxiliary gas. See Table 2 and Table 3 for additional information on using supplemental gas flows.

How Can I Introduce My Samples into the MS Detector?

You can introduce your samples into the MS detector in a variety of ways. Refer to Table 1.

The syringe pump is often used to introduce calibration solution for automatic tuning and calibrating in ESI mode. You can also use this technique to introduce a solution of pure analyte at a steady rate in ESI mode, for example, for determining the structure of an unknown compound.

You can also use a Tee union to direct samples from the syringe pump into an LC flow (without a column), which then enters the MS detector. This technique is used to introduce sample at a steady rate and at higher solvent flow rates; it is used especially for tuning in ESI or APCI on an analyte of interest. You can also use this technique to introduce a solution of pure analyte at a steady rate in ESI or APCI.

You can introduce samples from a syringe into the loop of the injector valve. You can then use the divert valve to introduce the sample into an LC flow, which then enters the MS detector. This technique is used in ESI or APCI to introduce pure analytes into the MS detector in a slug. It is useful when you have a limited quantity of pure analyte.

You can also use an LC autosampler to introduce samples into an LC flow. This technique is also used in ESI or APCI to introduce a slug of pure analyte into the LC flow and then into the MS detector.

Finally, you can perform LC/MS experiments by using an LC autosampler to introduce a mixture onto an LC column. This technique is used with ESI or APCI to separate the analytes before they are introduced sequentially into the MS detector.

You can refer to subsequent chapters in this manual and to *LTQ XL Getting Connected* for plumbing diagrams and methods of sample introduction.

	Sample Introduction Technique	Analytical Technique	Figure Reference
Syringe Pump Flow (no LC Flow)	Syringe pump*	ESI automatic tuning and calibrating ESI analysis of a pure analyte solution	<i>LTQ XL Getting Started</i> Figure 2-5

 Table 1.
 Sample introduction techniques

	Sample Introduction Technique	Analytical Technique	Figure Reference
LC Flow Without Chromatographic Separation (no column)	Syringe pump into LC flow (connected by Tee union)*	ESI or APCI automatic optimization of tuning on analyte of interest ESI or APCI analysis of a pure analyte solution	<i>LTO XL Getting Started</i> Figure 4-1 (ESI) Figure 6-1 (APCI)
	Loop injection into LC flow	ESI or APCI analysis of a pure analyte solution	<i>LTQ XL Getting Started</i> Figure 5-6 (ESI) Figure 8-1 (APCI)
	Autosampler injection into LC flow (one or multiple injections)	ESI or APCI analysis of a pure analyte solution	<i>LTQ XL Getting Connected</i> Figure 11-5 (ESI) Figure 11-8 (APCI)
LC Flow With Chromatographic Separation	Autosampler injections into LC column via LC flow (one or multiple injections	ESI or APCI analysis of mixtures	

Table 1.Sample introduction techniques

*Provides steady state introduction of sample (direct infusion)

What Types of Buffers Should I Use? What Types Should I Avoid?

Many LC applications use nonvolatile buffers such as phosphate and borate buffers. It is best to avoid the use of nonvolatile buffers with the MS detector because they can cause the following problems:

- Blocking the capillary in the probe
- Causing salt buildup on the spray head and thus compromising the integrity of the spray

Use volatile buffers when you use the MS detector. Many volatile buffer solutions are available that can be used instead of nonvolatile ones. Volatile buffer solutions can include the following:

- Acetic acid
- Ammonium acetate
- Ammonium formate
- Ammonium hydroxide
- Triethylamine (TEA)
- Trifluoroacetic acid

How Should I Set Up the MS Detector for Various LC Flow Rates?

The ESI probe can generate ions from liquid flows² of 1 μ L/min to 1.0 mL/min. This flow rate range allows you to use a wide range of separation techniques: CE, CEC, capillary LC, microbore LC, and analytical LC.

The APCI probe can generate ions from liquid flows³ of 200 μ L/min to 2.0 mL/min. This flow range allows you to use microbore LC, analytical LC, and semi-preparative LC.

As you change the rate of flow of solvents entering the MS detector, you need to adjust several of the MS detector parameters, as follows:

For ESI, you need to adjust the capillary temperature and adjust the gas flow rates for the Sheath, Auxiliary, and/or Sweep gas.

For APCI, you need to adjust the capillary temperature and vaporizer temperature and adjust the gas flow rates for the Sheath, Auxiliary, and/or Sweep gas.

In general, an increase in the rate of liquid flowing into the MS detector, requires a higher temperature of the ion transfer capillary (and vaporizer) and the higher gas flow rate.

Table 2 provides guidelines for ESI operation for ion transfer capillary temperatures and gas flow rates for various LC solvent flow rates.

Table 3 provides guidelines for APCI operation for the ion transfer capillary temperature, vaporizer temperature, and gas flow rate for a range of LC solvent flow rates.

² The ESI probe can generate ions from liquid flows of as low as 1 μ L/min. However, flows below 5 μ L/min require more care, especially with the position of the fused silica sample tube within the ESI probe.

 $^{^3}$ For the APCI probe, flows below 200 $\mu L/min$ require more care to maintain a stable spray.

1 Introduction

How Should I Set Up the MS Detector for Various LC Flow Rates?

LC Flow Rates	Suggeste d Column Size	lon Transfer Capillary Temperatu re	Sheath Gas	Auxiliary and/or Sweep Gas
Infusion or LC at flow rates of < 10 μL/min	Capillary	Typical setting: 150 to 200 °C	Not required Typical setting: 5 to 15 units	Not required Typical setting: 0 units
LC at flow rates from 50 to 200 µL/min	1 mm ID	Typical setting: 200 to 275 °C	Required Typical setting: 20 to 40 units	Not required, but might help depending on conditions Typical setting: 0 to 20 units
LC at flow rates from 100 to 500 µL/min	2 to 3 mm ID	Typical setting: 250 to 350 °C	Required Typical setting: 40 to 60 units	Not required, but usually helps to reduce solvent background ions Typical setting: 0 to 20 units
LC at flow rates from 0.4 to 1 mL/min	4.6 mm ID	Typical setting: 300 to 400 °C	Required Typical setting: 60 to 100 units	Required Typical setting: 10 to 40 units

5 1 51 , ,	Table 2.	Guidelines for setting operating parameters for LC/ESI/MS*
------------	----------	--

* Note: Be sure to choose either Auxiliary gas and/or Sweep gas according to the hints in Should I Use Sheath, Auxiliary, and/or Sweep Gases?

Table 3.	Guidelines for setting operating parameters for LC/APCI/MS*
----------	---

LC Flow Rate	lon Transfer Capillary Temperatur e	Vaporizer Temperatur e	Sheath Gas	Auxiliary and/or Sweep Gas
LC at flow rates from 0.2 to 2 mL/min	Typical setting: 150 to 225 °C	Typical setting: 400 to 550 °C	Required Typical setting: 40 to 100 units	Not required, but usually helps to reduce solvent background ions Typical setting: 0 to 20 units

* Note: Be sure to choose either Auxiliary gas and/or Sweep gas according to the hints in Should I Use Sheath, Auxiliary, and/or Sweep Gases?

What is Tuning and Calibration of the MS Detector All About?

To optimize the performance of data acquisition on the LTQ XL MS detector, tune and calibrate in four steps:

- In ESI mode, you infuse a calibration solution into the MS detector at a steady rate of 5 μ L/min for several minutes. In Tune Plus, you observe the signal at *m*/*z* 195, the mass-to-charge ratio of caffeine in the calibration solution. Then, while observing the signal at *m*/*z* 195, you adjust probe positions and gas flows to achieve the greatest signal strength while still maintaining a stable spray of ions into the MS detector.
- Once you have established a stable spray of ions into the MS detector, tune the MS detector. In this step, you use the automatic tuning procedure in Tune Plus to ensure that the transmission of ions into the MS detector is optimum. You observe the Tune Plus window as the Xcalibur data system tunes your LTQ XL MS detector automatically.
- After your tune method is optimized, calibrate the MS detector. In this step, you want to ensure that the calibration parameters complete automatic calibration successfully. The Calibrate dialog box in Tune Plus provides a readback of the status of the calibration parameters, both during the automatic calibration and when calibration is complete.
- Lastly, if you want to maximize the detection of one or more particular ions, you can optimize the tune of the MS detector with your analyte of interest in the ionization mode that you are going to use to analyze your samples. You choose a mass-to-charge ratio of your analyte of interest. Alternatively, you can choose an ion in the calibration solution that is closest to the mass-to-charge ratio for your ion of interest. (It is sometimes possible to acquire qualitative data without optimizing the parameters, but detection sensitivity might be compromised.)

Calibration parameters are instrument parameters whose values do not vary with the type of experiment. It is recommended that you calibrate the MS detector at least once every three months and that you check the calibration about once a week.

Automatic and semi-automatic calibration (including checking the calibration) require that you introduce calibration solution into the MS detector at a steady flow rate while the procedure is running. You introduce the solution directly from the syringe pump into the MS detector in the ESI/MS mode.

Tune parameters are instrument parameters whose values can vary with the type of experiment. For example, if your experiment requires quantitative data on one or more particular ions, you need to tune the MS detector with your analyte if you change any one of the parameters specific to the experiment or analyte.

Automatic and semi-automatic tuning procedures (including optimizing the collision energy) require that you introduce calibration solution, or a tuning solution of your analyte of interest, into the MS detector at a steady rate in either of two ways:

- Introduce the solution directly from the syringe pump. See Setting Up the Syringe Pump for Tuning and Calibration in Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode".
- Introduce the sample from the syringe pump into the effluent of the LC by using a Tee union. See Setting Up to Introduce Sample by Syringe Pump into Solvent Flow from an LC in Chapter 4: "Tuning with Your Analyte in LC/ESI/MS Mode".

Use the first method for tuning if you intend to use an experiment type at a low flow rate involving the syringe pump. The second method is useful if you intend to use an experiment type at a higher flow rate involving the LC. However, the second method of introduction puts a comparatively large amount of analyte into the MS detector. Therefore, before you can perform an analytical run to analyze for the analyte, you might need to clean the API spray shield.



CAUTION Do not use the calibration solution at flow rates above 10 μ L/min. Ultramark 1621 can contaminate your system at high concentrations.

In most cases, you can use the tune you obtain from the automatic or semi-automatic tuning procedures for your analytical experiments. However, for some applications, you might need to tune several MS detector parameters. In that case, you would tune manually. With the manual tuning process, you introduce a tuning solution at a steady flow rate.

Note The most important parameters that affect the signal quality during ESI/MS operation are the ion transfer capillary temperature, tube lens voltage, gases, and solution flow rate. For optimum sensitivity, tune with the instrument in the same operational mode as the mode you use for the analytical experiment.

Table 4 summarizes methods of sample introduction for each of the calibration and tuning procedures.

Calibrating			Tuning				
Sample/ Sample Intro	Check	Auto	Semi- auto	Auto	Semi- auto	Manual	Collision Energy
Calibration solution/ Syringe pump	✓	~	~	~	✓	✓	✓
Your tune solution/ Syringe pump				~	✓	~	✓
Your tune solution/ Syringe pump into LC flow by using Tee union				✓	~	✓	~

 Table 4.
 Summary of methods of sampleintroduction for calibration and tuning

What Types of Experiments Can I Perform with the LTQ XL MS Detector?

This topic describes several types of experiments that you can perform with the LTQ XL MS detector. The experiments can be grouped into the following categories:

- General MS or MSⁿ
- Data-Dependent[™]
- Ion Mapping[™]
- Ion Tree

You can specify which type of experiment you want to perform in the Instrument Setup window, and then save it in an Instrument Method (.METH) file.

Note Procedures for these experiments are beyond the scope of this *LTQ XL Getting Started* manual. If you need more information, refer to online Help.

You can use a General experiment to collect qualitative data for structural analysis. The Xcalibur data system includes an Instrument Method template in Instrument Setup so you can get started with a General MS or MSⁿ experiment. For an example of a General MS or MSⁿ experiment template, see Figure 1

In a General *MS* experiment, you need to specify the mass range of your analyte(s) of interest. In a General *MS/MS* experiment, you need to specify a parent (precursor ion) that fragments into distinctive product ions. In a General MS^n experiment, you need to specify the mass-to-charge ratios of *all* the parent ions of interest. The LTQ XL MS detector can then collect data on the ions in the range or on the product ions of the parent ion(s) that you specify.

If you use a General experiment to collect data for qualitative (structural) analysis, you specify the scan mode (MS through MSⁿ) for which you want data in the Scan Event Settings group box. If you specify MS/MS or MSⁿ, you then choose the parent ion(s) for which you want data in the MSⁿ Settings table. The LTQ XL MS detector can then collect distinct qualitative information for structural analysis or for spectral reference.

General MS or MSn Experiments

General MS or MSn



Figure 1. MS Detector Setup page in Instrument Setup, showing a template for a General MS experiment

The LTQ XL MS detector can generate reproducible, analyte-specific spectra, even from laboratory to laboratory. Consequently, reference spectra that are generated with the LTQ XL MS detector can be used to confirm structures of compounds generated with other LTQ XL systems.

Data-Dependent Experiments

Data dependent <u>M</u>S/MS

Data dependent triple play

A Data-Dependent experiment is best used for the qualitative analysis of unknown compounds for structure elucidation or confirmation. The LTQ XL MS detector uses the information in a data-dependent experiment to make decisions about the next step of the experiment automatically—without input from a user. Instrument Setup contains the Instrument Method templates that you need to get started with data-dependent experiments. For an example of a data-dependent Triple Play experiment template, see Figure 2. A data-dependent experiment produces a great deal of data from a single sample analysis. You can run a data-dependent experiment even if you know very little about your sample, and even if you are unfamiliar with the variables of mass spectroscopy. In a data-dependent experiment, you can specify parent ions for fragmentation or you can let the LTQ XL MS detector automatically select the ions for fragmentation. The LTQ XL MS detector can collect the structural information for every parent ion in the sample automatically, even if the sample is a mixture of compounds.

A data-dependent experiment requires minimal input from a user about how the experiment should best proceed. The user specifies that one or more scan events of an experiment segment are to be run as data-dependent. Then, the LTQ XL MS detector collects MS/MS or MSⁿ data and makes decisions about what the next step in the experiment should be to collect even more data. For example, in a data-dependent Triple Play experiment for a mixture of compounds, the LTQ XL MS detector can decide which parent ion to isolate, the charge state of the parent ion, and the molecular mass of the compound.

Ion Mapping experiments can be data-dependent. (The Total Ion Map, Neutral Loss Ion Map, and Parent Ion Map experiments are *not* data-dependent.) The Data-Dependent Zoom Map experiment collects ZoomScan data on every scan interval in a specified mass range.

Ion Tree experiments are types of data-dependent experiments. These experiments provide methods for automatically interpreting MSn data and arranging the data in formats that are easy to manipulate.

You can approach the setup of data-dependent experiments in either of two ways:

• If you have some idea of the parent ion, or if you expect a certain kind of parent, you can set up a list of possible parent ions. Then, when one of the parent ions you specified is detected, you can acquire product spectra and analyze the information. Conversely, you can also set up a list of ions that you do not want to be selected for fragmentation.



- **Figure 2.** MS Detector Setup page in Instrument Setup, showing a template for a Data-Dependent Triple Play experiment. (To select a scan event that makes active the Dependent Scan check box, click either the Scan Event 2 or Scan Event 3 button.
 - If you have little information about your compound, you can set up the parameters of a data-dependent experiment so that if the intensity of the ion signal is above a specified threshold, the LTQ XL MS detector generates product spectra. Parameters that you might specify, for example, include threshold values for the intensity of the MS or MSⁿ ion signal. Whatever threshold values you choose should accomplish the isolation of your parent ions of interest.

You can find useful structural information about your compound automatically with the simplest data-dependent experiment, Data-Dependent MS/MS. You specify the MS scan range, and you do not even need to specify a parent ion. The LTQ XL MS detector can then collect full scan MS data, pick the most intense parent ion in the spectrum, and fragment the ion to generate product ions.

A Data-Dependent Triple-Play experiment is the same as Data-Dependent MS/MS, but includes the identification of the charge state of the parent with the LTQ XL ZoomScan feature. A Data-Dependent Triple-Play experiment collects full scan MS data, and then uses ZoomScan to determine the charge state of the parent ion and calculate the molecular mass. The parent ion is then fragmented into product ions (MS/MS). For example, if the LTQ XL MS detector determines a charge state equal to 2, and if the mass-to-charge ratio of the parent ion is m/z 500, then the mass-to-charge ratios of the product ions can be up to m/z 1000 (or 2 × 500).

Use a data-dependent experiment (from templates in Instrument Setup) to do the following:

- Identify low-level impurities in high-purity compounds (Data-Dependent MS/MS)
- Identify metabolites in a complex mixture (Chromatographic Separation with Data-Dependent MS/MS)
- Build a custom library of composite MSⁿ spectra (Ion Tree)

You can use a Data-Dependent MSⁿ experiment to identify process impurities. In the quality assurance process for aspirin, for example, the LTQ XL MS detector can identify impurities of less than 0.1%.

A Data-Dependent MS/MS experiment of a complex mixture of drug metabolites can provide highly specific structural information. Characteristic masses along the metabolic pathways of a drug, for example, can produce MS/MS spectra that are specific to the structure of the drug. These spectra are essential in metabolite identification.

A data-dependent experiment can produce a composite spectrum of, for example, MS2, MS3, and MS4 data. The LTQ XL MS detector can store the MSⁿ fingerprint data in a custom MSⁿ library spectrum. The data is valuable for use in process control, quality assurance, or research.

Ion Mapping Experiments

Ion mapping...

An Ion Mapping experiment is best used to get full structural characterization of unknown molecules in complex mixtures. In an Ion Mapping experiment, you can get product ion scans on every parent ion over a specified mass range. An Ion Mapping experiment can help to identify automatically which parent ions were fragmented to yield a specified product ion. The experiment "maps" one or more parent ions by using the information from product ion scans.

The LTQ XL MS detector includes the following Ion Mapping templates in Instrument Setup so you can get started with an Ion Mapping experiment:

- Total (or full scan) Ion Map
- Neutral Loss Ion Map
- Parent Ion Map

These Ion Mapping experiments, in general, require that sample solution enter the MS Detector at a composition that is constant throughout. Therefore, you use infusion to introduce your sample for these Ion Mapping experiments. See Figure 3 for an example of an Ion Mapping experiment template.

Select ion map type	1
T <u>o</u> tal ion map	
Neutral Joss ion map	
Parent ion map	
	1

1 Introduction

What Types of Experiments Can I Perform with the LTQ XL MS Detector?

Untitled - Instrumen	nt Setup		_O×
Eile LTQXL Help	?		
Aglent1100 Bin	Total Ion Map Syringe Pump Divert Valve Contact Closure Summ Acquire time (min): Mass range:	100.00 Normal Positive C:Wcalibur\methods\Default_ESI.LTQT 300.00-2000.00	
'LTQ XL MS Surveyor LC Pump Thermo Pal	Parent mass gtep size (m/z): [solation width (m/z): Normalized goliision energy (%): Activation Q: Activation time (msec): Product mass rgnge (m/z):	1.0 2.0 36.0 0.250 30.000	
Ready	New method	Tyne Plus Help	NOT SAVED

Figure 3. Total lon Map page in Instrument Setup, showing a template that contains parameters for an Ion Mapping experiment

In a Total (or full scan) Ion Mapping experiment, you get product ion scans for each parent ion, so you can determine which parent ions lost a particular fragment to yield a particular product ion. Furthermore, you can determine which parent ions are related to specific product ions. For example, you can map the spectral peaks in a mass range from m/z 400 to m/z 2000 and specify to scan for MS/MS product ions in incremental steps of every mass-to-charge ratio, every fifth mass-to-charge ratio, or every tenth mass-to-charge ratio.

A Neutral Loss Ion Mapping Experiment collects scans for masses that have lost neutral fragments. As with Full Scan Ion Mapping, you can get product ion scans on every parent ion. However, a Neutral Loss Ion Map identifies which parent ions lost a neutral fragment of a particular mass. For example, you can specify a neutral loss of 80 u (as in the case of a phosphorylated peptide in a tryptic digest). A Neutral Loss Ion Mapping experiment can step through each product mass in the mixture. The experiment searches for evidence of the loss of a neutral moiety of mass 80 u.

A Parent Ion Mapping experiment identifies all the ions that produce a particular molecular ion that you specify. For example, if you specify a product ion mass of m/z 50, a Parent Ion Map includes all the parent ions that yielded the specified product ion, m/z 50.

A Data-Dependent Zoom Map is an Ion Mapping experiment that collects ZoomScan data on every scan interval in a mass range that you specify, as well as Data-Dependent MS/MS product spectra on every mass above an intensity threshold.

The results of any of the Ion Mapping experiments can be viewed in the Xcalibur Qual Browser window.

Ion Tree Experiments

Data dependent ion tree

In an Ion Tree experiment, the LTQ XL MS detector can collect MSⁿ data automatically. You can specify a particular parent ion for fragmentation, or you can let the LTQ XL MS detector find the parent ions automatically and fragment them to any level between MS² and MS¹⁰. The LTQ XL MS detector automates the collection of data by deciding what actions need to occur next for the experiment to progress. See Figure 4 for an example of an Ion Tree experiment template.

1 Introduction

What Types of Experiments Can I Perform with the LTQ XL MS Detector?

File IT 024. beb Image: It 024. beb Data Dependent Ion Tree Springe Pump Directivated Closure Summary Image: It 024. beb Data Dependent Ion Tree Springe Pump Directivated Closure Summary Image: It 024. beb Data Dependent Ion Tree Springe Pump Directivated Closure Summary Image: It 024. beb Image: Itom Closure Image: Itom C	gonacieu - inscrumer	it setup	
Image: Source of the set	<u>File L</u> TQ XL <u>H</u> elp		
Data Dependent Ion Tree Syinge Punc. Divent Valve Contact Closure Summary Agilent 100 Bin Experiment setting: Excur: © Breadh © Depth Max gendent: 5 Max gendent: 5 Analysis mode: © Infusion © Chromotographic Fun setting: Acquire time (min): 100.00 Turg method CVcsatou/methods/Default_ESILITOT		8	
New method Tune Help	Agilent1100 Bin Agilent1100 Bin "LTQ XL MS Surveyor LC Pump Thermo Pal	Data Dependent Ion Tree Syringe Pane Direct Valve Contact Closure Summary Experiment settings Ecour: © Breadth Deph Max geptit: 3	
	Ready	NOT SAVED	1

Figure 4. Data-Dependent Ion Tree page in Instrument Setup, showing a template for an Ion Tree experiment

In an Ion Tree experiment, you can specify either of two options that prioritize how the LTQ XL MS detector gathers information: depth focus and breadth focus.

- Depth focus characterizes an ion by performing a series of MSⁿ-level fragmentations (for example, MS/MS, MS3, MS4, etc.) before characterizing the next most intense ion in the MSⁿ series.
- Breadth focus characterizes all ions to the same MSⁿ level before advancing to the next MSⁿ level.

For example, if you specify a Maximum Depth of 3 and a Maximum Breadth of 2 in an Ion Tree experiment, the following occurs.
First, with either depth or breadth focus, the LTQ XL MS detector scans for parent ions (MS) over the specified mass range. The most intense ion of the MS spectrum is selected for fragmentation (MS/MS).

- Second, if you chose the depth focus, after the most intense ion of the MS spectrum is fragmented—producing an MS/MS spectrum—the LTQ XL MS detector selects and fragments the most intense ion of the *MS/MS* spectrum. This results in an MS3 spectrum, the level specified as the maximum depth for this example. The LTQ XL MS detector then backs up one level and fragments the second most intense ion of the *MS/MS* spectrum, creating more product ions on the level of MS3 from this parent ion. This process is then repeated for the second most intense ion in the *MS* spectrum.
- If you chose the breadth focus, after the most intense ion of the MS spectrum is fragmented—producing an MS/MS spectrum—the LTQ XL MS detector selects and fragments the second-most intense ion of the *same* MS spectrum. The fragmentation of parent ions continues to the *Max Breadth* level that you specified (2, for this example). After the two most intense peaks on the MS level are fragmented, the LTQ XL MS detector scans the first *MS/MS* spectrum to select and fragment the two most intense ions. This results in product ions on the level of MS3, the level specified as the maximum depth for this example. This process is then repeated for the second most intense ion in the *MS* spectrum.

The results of a Data-Dependent Ion Tree experiment can be viewed in the Xcalibur Qual Browser window. The results are displayed as a structure tree that originates from a particular parent ion.

Chapter 2 Setting Up the Ion Source for Tuning and Calibrating the MS Detector

This chapter provides information on setting up the hardware for tuning and calibrating your LTQ XL MS detector. You tune and calibrate the MS detector in the ESI mode before you acquire data in either the ESI or APCI mode.

This chapter contains the following topics:

- Placing the LC/MS System in Standby
- Removing the APCI Probe
- Removing the Ion Max Ion Source Housing (optional)
- Installing the Ion Sweep Cone (optional)
- Installing the Ion Max Ion Source Housing
- Installing the ESI Probe

Placing the LC/MS System in Standby

	1	Programs 🕨		Accessor	ies⊧ ⊧	Ľ	
nal		<u>D</u> ocuments►		Xcalibur	,	8	Instrument Configuration
ssio	V -	Settings	Γ			•	LTQTune
rofe		Search 🕨				*	Merlin Xcalibur
XP P	•	Help				170	
dows		<u>R</u> un					
Win	0	Shut Down					
<u>a</u>	Start						





The LC/MS system needs to be placed in Standby condition before you can remove the ion source.

Place the LC/MS system in Standby

- 1. If necessary, stop the flow of solvent to the API source as follows:
 - a. If the Xcalibur data system is not already open, choose **Start** > **Programs** > **Xcalibur** > **Xcalibur** from the Windows[®] taskbar to open the Xcalibur window.
 - b. In the Xcalibur Home Page window Roadmap view, choose
 GoTo > Instrument Setup to open the Instrument Setup window.
 - c. Click the **Surveyor**[®] **MS Pump** button on the view bar in the Instrument Setup window to display the Surveyor MS Pump view.
 - d. Choose **Surveyor MS Pump > Direct Control** to open the Surveyor MS Pump Direct Control dialog box.
 - e. In the Direct Control dialog box, click the **Pump Off** button to stop the MS pump.
- 2. If Tune Plus is not already open, choose **Start > Programs > Xcalibur > LTQ XLTune** from the taskbar to open Tune Plus.

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

3. If the MS detector is On, click the **On/Standby** button to place the MS detector in the Standby mode. When the MS detector is in Standby, the LTQ XL MS detector turns off the ion source sheath gas, auxiliary gas, and high voltage.

The LC/MS system is now in Standby and it is safe to remove the ion source.

If the ESI probe is already installed in the Ion Max[™] ion source housing, leave the LC/MS system in Standby and go to Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode".

If the ESI probe is not already installed in the Ion Max ion source housing, go to the next section, Removing the APCI Probe.

Removing the APCI Probe

This topic describes how to remove the APCI probe from the Ion Max ion source housing.

Note The following procedures assume that you are familiar with your instrument and software. If you need additional guidance, refer to LTQ XL online Help, *LTQ XL Getting Connected, Ion Max API Source Hardware Manual*, or the *LTQ XL Hardware Manual*.



CAUTION AVOID BURNS. At operating temperatures, the APCI vaporizer can severely burn you! The APCI vaporizer typically operates between 400 and 600 °C. Always allow the heated vaporizer to cool to room temperature (for approximately 20 min) before you touch or remove this component.

Remove the APCI probe

- 1. Unplug the vaporizer heater cable from the vaporizer heater cable socket on the APCI probe. See Figure 5.
- 2. Disconnect the sample transfer line from the APCI probe. (See Figure 5.)
- 3. Remove the auxiliary gas line (green-colored fitting) from the APCI probe. (Figure 5)
- 4. Remove the sheath gas line (blue-colored fitting) from the APCI probe.



Figure 5. Ion Max ion source housing with APCI probe installed



CAUTION AVOID BURNS. At operating temperatures, the APCI vaporizer can severely burn you! The APCI vaporizer typically operates between 400 and 600 °C. Always allow the heated vaporizer to cool to room temperature (for approximately 20 min) before you touch or remove this component.

- 5. Remove the APCI probe as follows:
 - a. Connect the vaporizer heater cable to the ESI interlock socket on the ion source housing. See Figure 6.
 - b. Release the probe locking lever to loosen the probe collar. You might need to unscrew the lever a few turns to permit probe movement.
 - c. Carefully pull the probe straight back in the port in the housing until it meets with the slot in the API interlock block. The guide pin on the probe manifold will prevent you from rotating the probe until the pin is aligned with the slot in the API interlock block. Once the probe is all the way back and aligned with the slot, turn the probe 45 degrees counter-clockwise to free the probe from the alignment notch.
 - d. Pull the probe straight out to remove it from the ion source housing.
 - e. Store the APCI probe in its original shipping container.



Figure 6. Ion Max ion source housing, detail of components

- 6. Remove the 8 kV cable from the corona needle high voltage receptacle as follows:
 - a. Unlock the cable by rotating the locking ring counter-clockwise.
 - b. Unplug the 8 kV cable from the corona needle high voltage receptacle.



CAUTION AVOID INJURY. The corona discharge needle is very sharp and can puncture your skin. Handle it with care.

- 7. Remove the corona needle as follows:
 - a. Unlock the ion source housing door by turning the locks 90 degrees so that the knobs are horizontal.
 - b. Open the ion source housing door.
 - c. Using pliers, grasp the corona needle and pull it straight out of the corona needle contact. See Figure 7.

- d. Close and lock the ion source housing door.
- 8. Store the corona needle in its original shipping container.

The APCI probe and the corona needle are now properly removed from the Ion Max ion source housing.

If you want to install the optional ion sweep cone, go to the next section, Removing the Ion Max Ion Source Housing.

If you do not want to install the ion sweep cone, go to "Installing the ESI Probe" on page 39.



Figure 7. Corona needle, view from rear

Removing the Ion Max Ion Source Housing

The Ion Max ion source housing is removed to access the ion sweep cone.

Note If an ion source probe is still installed in the ion source housing, the external liquid lines should first be disconnected before removing the ion source housing.

Remove the ion source housing

- 1. Remove the drain tube from the ion source housing drain. See Figure 8.
- 2. Rotate the ion source housing locking levers 90 degrees to release the ion source housing from the ion source mount assembly.
- 3. Remove the ion source housing by pulling the housing straight off of the ion source mount assembly
- 4. Place the ion source housing in a safe location for temporary storage.

The Ion Max ion source housing is now properly removed.





Installing the lon Sweep Cone

The ion sweep cone is a metallic cone that is installed over the ion transfer tube. The ion sweep cone channels the sweep gas towards the entrance of the capillary. This helps to keep the entrance of the ion transfer tube free of contaminants. The net result is a significant increase in the number of samples that can be analyzed without a loss of signal intensity. In addition, keeping the ion transfer tube entrance cleaner reduces the need for frequent MS detector maintenance.

Install the ion sweep cone

- 1. Remove the ion sweep cone from its storage container. Inspect and clean it if necessary.
- 2. Note the location of the sweep gas supply port in the API cone seal. The gas inlet on the ion sweep cone is placed in this port. See Figure 9 and Figure 10.



Figure 9. Sweep gas supply port in the API cone seal



Figure 10. Ion sweep cone, showing the gas inlet



CAUTION AVOID BURNS. At operating temperatures, the ion transfer tube can severely burn you! The ion transfer tube typically operates between 200 and 400 °C. Always allow the ion transfer capillary to cool to room temperature (for approximately 20 min) before you install the ion sweep cone. Always be careful not to touch the entrance end of the ion transfer tube when it is exposed.

- 3. After the ion transfer tube has cooled to room temperature, carefully align the gas inlet on the ion sweep cone with the sweep gas supply port in the API cone seal. Firmly press the ion sweep cone into position.
- 4. If necessary to achieve a proper ion sweep cone installation, you might adjust the set screws around the perimeter of the ion sweep cone.

The ion sweep cone is now properly installed on the MS detector.

Installing the Ion Max Ion Source Housing

Reinstall the Ion Max ion source housing

1. Carefully align the two guide pin holes on the rear of the ion source housing with the ion source housing guide pins on the MS detector, and carefully press the ion source housing onto the ion source mount. See Figure 11 and Figure 12.



Figure 11. Rear view of the Ion Max ion source housing

2 Setting Up the Ion Source for Tuning and Calibrating the MS Detector Installing the Ion Max Ion Source Housing



Ion Source Housing Guide Pins



2. Rotate the ion source housing locking levers 90 degrees to lock the ion source housing onto the ion source mount assembly.



CAUTION Prevent solvent waste from backing up into the ion source and MS detector. Always ensure that liquid in the drain tube is able to drain to a waste container.

3. Reinstall the ion source drain tube as follows:



CAUTION Do **not** vent the API source drain tube (or any vent tubing connected to the waste container) to the same fume exhaust system to which you have connected the forepumps. The analyzer optics can become contaminated if the API source drain tube and the (blue) forepump exhaust tubing are connected to the same fume exhaust system.

CAUTION Your laboratory must be equipped with at least two fume exhaust systems. Route the (blue) forepump exhaust tubing to a dedicated fume exhaust system. Route the drain tube from the API source to a waste container. Vent the waste container to a dedicated fume exhaust system.

- a. Connect the 1-in. ID Tygon' tubing to the ion source housing drain.
- b. Attach the free end of the hose to a dedicated drain system. Ideally, the drain system should be vented to a fume exhaust system.

The Ion Max ion source housing is now properly installed on the MS detector.

Installing the ESI Probe

Install the ESI probe

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

Note If your ESI probe does not already have a sample tube (fused-silica capillary or metal needle) and safety sleeve attached, follow the procedure for installing a sample tube and PEEK safety sleeve that is outlined in **Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve** in the *Ion Max API Source Hardware Manual*.

- 2. Ensure that the probe locking lever on the ion source housing is unlocked (opened to its widest position). See Figure 13.
- 3. Insert the ESI probe into the port in the ion source housing, align the guide pin on the probe body at a minus 45 degree angle from the API interlock block. See Figure 14



Figure 13. Ion Max ion source housing probe locking lever open



Figure 14. ESI probe, side view

- 4. Push the probe into the port until the guide pin meets with the probe collar on the ion source housing.
- 5. Turn the probe 45 degrees clockwise and align the guide pin with the slot in the API interlock block (you might need 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 rotating the probe locking lever towards the front of the housing; closing the probe locking lever towards the rear of the ion source housing might make it difficult to unlock. You might first need to tighten the locking lever threaded shaft by rotating it clockwise a few turns if rotating the lever does not tighten the probe collar enough.
- 7. Insert the APCI vaporizer heater cable into the API interlock socket.



8. Insert the stainless steel ZDV fitting (grounding union) into the grounding bar on the ion source housing. See Figure 15.

Figure 15. Ion Max ion source housing with ESI probe installed

- 9. Connect the sheath gas fitting (blue) to the sheath gas inlet (S) on the probe. (See Figure 15.)
- 10. Connect the auxiliary gas fitting (green) to the auxiliary gas inlet (A) on the probe. (See Figure 15.)
- 11. Connect the 8 kV cable to the ESI needle high voltage receptacle on the ESI probe. Tighten the locking ring on the 8 kV connector.
- 12. Connect the sample transfer tubing to the grounding union.

The ESI probe is now properly installed in the Ion Max ion source housing.

Leave the LC/MS system in Standby and go to Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode".

Chapter 3 Tuning and Calibrating Automatically in the ESI/MS Mode

This chapter provides information on how to tune and calibrate the LTQ XL MS detector in the ESI/MS mode. For most applications, you tune and calibrate in the ESI mode through automatic procedures. The procedures use a calibration solution that is introduced into the MS detector in low flow mode. The procedures properly tune and calibrate the MS detector for ESI operation (refer to Table 2 on page 2 for ESI operating parameter guidelines). You need to calibrate the MS detector every one to three months of operation for optimum performance over the entire mass range of the detector.

To tune and calibrate your MS detector automatically in the ESI/MS mode, you do the following:

- Infuse a low concentration calibration solution containing caffeine, MRFA, and Ultramark 1621 into the ESI source by using the syringe pump. (Refer to the section, Setting Up the Syringe Pump for Tuning and Calibration.)
- Test the efficiency and stability of the spray of calibration solution into the MS detector. You can observe the following singly-charged, positive ions for caffeine, MRFA, and Ultramark 1621 in the Tune Plus window: *m*/*z* 195, 524, 1222, 1522, and 1822.
- Tune the MS detector from the Tune Plus window to optimize automatically the lenses.
- Calibrate the MS detector to adjust automatically the voltages of the linear trap.

This chapter contains the following sections:

- Setting Up the Syringe Pump for Tuning and Calibration
- Setting Up the MS Detector in the Xcalibur Data System for Tuning and Calibration
- Testing the Operation of the MS Detector in the ESI/MS Mode

- Tuning the MS Detector Automatically in the ESI/MS Mode
- Saving Your ESI/MS Tune Method
- Calibrating the MS Detector Automatically
- Cleaning the MS Detector after Tuning and Calibrating

Setting Up the Syringe Pump for Tuning and Calibration

You introduce tuning and calibration solution into the API source with a syringe infusion pump. A syringe pump allows you to infuse a sample solution into the API source for extended periods of time.

The syringe pump and syringe are located on the front panel of your LTQ XL MS detector. To infuse solution for tuning and calibration, you install on the pump a 500-mL Unimetrics[®] syringe containing the calibration solution.

Note To minimize the possibility of cross-contamination, use a different syringe and section of fused silica tubing for the calibration solution than you do for your sample solution.

Set up the syringe pump for infusion

1. Connect a 4 cm (1.5 in.) segment of Teflon[®] tube with a (brown) fingertight fitting and a (brown) ferrule to the (black) LC union. See Figure 16.



Figure 16. Plumbing connections for the syringe

- 2. Load a clean, 500-μL Unimetrics syringe with 450 μL of the calibration solution. (Refer to Appendix A: "Sample Formulations" on page 139 for a procedure for making the calibration solution.)
- 3. Insert the syringe needle into the segment of Teflon tube.
- 4. Place the syringe into the syringe holder of the syringe pump.
- 5. While squeezing the blue release button on the syringe pump handle, push the handle forward until it just contacts the syringe plunger.
- 6. Connect a fused-silica infusion line from the LC union to the (stainless steel) grounding union as follows. See Figure 17.

- a. Connect the infusion line with a (brown) fingertight fitting and a (brown) ferrule to the free end of the LC union.
- b. Connect the other end of the infusion line with a (red) fingertight fitting and a (brown) ferrule to the grounding union.

The syringe pump is now properly set up for infusing solution into the MS detector.

Go to the next section, Setting Up the MS Detector in the Xcalibur Data System for Tuning and Calibration.



Figure 17. ESI/MS plumbing connections for the fused-silica infusion line

Setting Up the MS Detector in the Xcalibur Data System for Tuning and Calibration



You first tune manually with calibration solution to establish a stable spray of solution and to ensure that enough ions are detected to calibrate the MS detector. You then calibrate the MS detector automatically to optimize the parameters that affect ion detection. With the optimized MS detector, the Xcalibur data system can isolate and fragment ions and determine their mass-to-charge ratios. Perform a calibration periodically, every one to three months, for optimum performance of the MS detector.

Note The following procedures assume that you are familiar with your LTQ XL instrument and the Tune Plus window. If you need additional guidance, see LTQ XL online Help, *LTQ XL Getting Connected*, and/or the *LTQ XL Hardware Manual*.

CAUTION Before you begin normal operation each day, ensure that you have sufficient nitrogen for your API source. If you run out of nitrogen, the LTQ XL MS detector automatically turns Off to prevent the possibility of atmospheric oxygen from entering the ion source. The presence of oxygen in the ion source when the MS detector is On could be unsafe. (In addition, if the LTQ XL MS detector automatically turns Off during an analytical run, you could lose data.)

Set up the MS detector in the Xcalibur data system for tuning and calibration in the ESI/MS mode

- 1. If you have not already done so, open the Tune Plus window from the Start button on your Windows XP task bar, as follows:
 - a. Choose **Start > Programs > Xcalibur > Xcalibur** to display the Xcalibur Home Page Roadmap view.
 - b. Click the **Instrument Setup** button to display the Instrument Setup window.
 - c. Click the **Finnigan LTQ XL** button to display the New Method page.
 - d. Click the **Tune Plus** button to display the Tune Plus window. See Figure 18.
- 2. In the Tune Plus window, on the Control/Scan Mode toolbar, click the **On/Off/Standby** button to take the MS detector out of the Standby (or





Setting Up the MS Detector in the Xcalibur Data System for Tuning and Calibration

Off) mode and turn it On. When you turn the MS detector to On, you initiate the following events:

- The MS detector begins scanning.
- Nitrogen flows into the ESI probe.
- The LTQ XL MS detector applies high voltage to the ESI probe.
- The Xcalibur data system shows a real-time display in the Spectrum view.



Figure 18. Tune Plus window, showing the MS detector in the Standby mode

Note The Xcalibur data system contains customized tune files for different applications in the folder *C:\Xcalibur\methods*, including one for low flow LC/ESI/MS operation.

- 3. Open the Tune Method file that stores the factory default tune settings for low-flow ESI operation, as follows:
 - a. Choose **File > Open** to display the Open dialog box.
 - b. Browse for the folder C:\Xcalibur\methods. Select the file AutoTune.LTQTune.
 - c. Click **Open** to open the file. Tune Plus downloads the Tune Method parameters to the MS detector.
- 4. Examine the pre-tune ESI source settings as follows:
 - a. From the Instrument Setup toolbar, click the **API Source** button to open the ESI Source dialog box. Verify that the settings in your dialog box are the same as those shown in Figure 19.
 - b. Click **OK** to return to the Tune Plus window.

ESI Source	×
	Actual
Sheath Gas Flow Rate (arb): 10	9.90
Aux <u>G</u> as Flow Rate (arb): 0	0.01
S <u>w</u> eep Gas Flow Rate (arb): 0	0.00
Spray Voltage (kV) : 4.50	4.50
Spray Current (µA):	0.05
Capillary Temp (*C): 275.00	274.99
Capillary Voltage (V): 13.00	13.04
Tube Lens Offset (V): 100.00	100.15
Apply OK Cancel <u>H</u>	elp

- Figure 19. ESI Source dialog box, showing the settings to start a typical low flow experiment
- 5. Set the scan parameters for tuning and calibration, as follows:



3 Tuning and Calibrating Automatically in the ESI/MS Mode

Setting Up the MS Detector in the Xcalibur Data System for Tuning and Calibration



- a. On the Control/Scan Mode toolbar, click the **Define Scan** button to open the Define Scan dialog box. See Figure 20. (If your dialog box appears different from the one shown in the figure, it is probably because the advanced settings are not displayed. You can turn on the advanced settings as follows: In Tune Plus, choose **ScanMode**, and then click **Advanced Scan Features** to select the option.)
- b. In the Scan Description group box, in the Mass Range list box, select Normal to allow for a selection of mass ranges between m/z 150 to 2000.
- c. In the Scan Rate list box, select Normal to specify a normal scan rate.
- d. In the Scan Type list box, select Full specify a full scan.
- e. In the Scan Time group box, in the Microscans spin box, enter 1 to set the total number of microscans to 1.
- f. In the Max. Inject Time spin box, enter 200.000 to specify a 200 ms maximum injection time.
- g. In the Source Fragmentation group box, confirm that the On check box is not selected (
) to specify that the ion source fragmentation option is turned off.
- h. In the Scan Ranges group box, in the Input list box, select From/To to make available the First Mass and Last Mass text boxes in the Scan Ranges table.

Setting Up the MS Detector in the Xcalibur Data System for Tuning and Calibration

Define Scan											×
Scan History: ITMS + c Full ms [150.00-2000.00]											
Scan Description	n Description MSn Settings						Scan Ranges				
<u>M</u> ass Range: Normal ▼ Scan Bate: Normal ▼	n	Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)		#	First Mass (m/z)	Last Mass (m/z)	
	2		1.0	20.0	0.250	10.000		1	150.00	2000.00	
Scan Lype: Full	3		1.0	20.0	0.250	10.000		2			
	4		1.0	20.0	0.250	10.000		3			
	5		1.0	20.0	0.250	10.000		4			
Scan Time	6		1.0	20.0	0.250	10.000		5			
Mi <u>c</u> roscans: 1	7		1.0	20.0	0.250	10.000		6			
Max Inject Time (ms): 200,000	8		1.0	20.0	0.250	10.000		7			
	9		1.0	20.0	0.250	10.000		8			
	10		1.0	20.0	0.250	10.000		9			
Source Fragmentation							10 _!	put: From/	To 💌		
	Арр	ly	OK	Cancel	Hel	р	In	iectic	n B <u>E</u>	Acti <u>v</u> ation]

Figure 20. Define Scan dialog box, showing the default settings for ESI/MS operation

- i. In the Scan Ranges group box, in the Scan Ranges table, in the First Mass text box, enter 150 to set the first mass for the scan range to m/z 150.
- j. In the Last Mass text box, enter 2000 to set the last mass for the scan range to m/z 2000.
- k. Ensure that the settings in your Define Scan dialog box are the same as those shown in Figure 20.
- 1. Click **OK** to apply the MS detector scan parameters and to close the Define Scan dialog box.
- 6. On the Control/Scan Mode toolbar, click the **Centroid/Profile** button to toggle the data type to profile. (The picture on the button should be the same as that shown here.)
- 7. Click the **Positive/Negative** button to toggle the ion polarity mode to positive. (The picture on the button should be the same as that shown here).

The MS detector is now properly set up in the Xcalibur data system for tuning and calibration in the ESI/MS mode.





Testing the Operation of the MS Detector in the ESI/MS Mode

You are now ready to test whether your MS detector is operating properly. To test for proper operation, you infuse the calibration solution into the ESI source, and then you monitor the real-time display of the mass spectrum of calibration solution. You want to ensure that a stable spray of solution enters the MS detector.

Test the operation of the MS detector in the ESI/MS mode



1. Click the **Syringe Pump** button to display the Syringe Pump dialog box. See Figure 21.

Syringe Pump				×
Flow Control				
				Actual
⊙ <u>O</u> n	Flow \underline{R} ate (µL/min):	5.00	÷	0.00
O O <u>f</u> f				
Туре				
C Hamilton	Volume (uL):	500	-	
<u> <u> unimetrics</u> </u>	Suringe ID (mm):	3.260		
C Other	olimiliao Tor (uuu):	1	<u>~</u>	
·				
Apply	UK	Cancel		Help

Figure 21. Syringe Pump dialog box

- 2. Turn on the syringe pump and set an infusion flow rate of 5 $\mu L/\text{min},$ as follows:
 - a. In the Flow Control group box, click the **On** option button to make active the Flow Rate spin box.
 - b. Type 5 in the Flow Rate spin box to specify a rate of 5 μ L/min.

Note This procedure assumes that you are using the $500-\mu L$ Unimetrics syringe that is provided with your LTQ XL system. If you are using another type of syringe, select the option button corresponding to your syringe.

c. If you are using a standard Unimetrics (or Hamilton) syringe, set up the syringe parameters as follows:

- i. In the Type group box, click the **Unimetrics** (or **Hamilton**) option button to specify the proper syringe type.
- ii. Click the **Volume** list box arrow to display the list of available volumes, and then select *500* (or your syringe size) from the list to set the proper syringe volume. Note that, if you are using a Unimetrics syringe, the LTQ XL MS detector automatically sets the syringe ID to its proper value of 3.260 mm.
- d. If you are not using a Unimetrics (or Hamilton) syringe, set up the syringe parameters as follows:
 - i. In the Type group box, click the **Other** option button to make active the syringe ID spin box.
 - ii. Type the inner diameter of your syringe in the Syringe ID spin box.
- e. Click **OK** to apply the syringe parameters, start the syringe pump, and close the Syringe Pump dialog box.
- 3. On the File/Display toolbar, click the **Display Spectrum View** button to ensure that the Spectrum view is displayed.
- 4. Monitor the data for the calibration solution, as follows:
 - a. In the Spectrum view of the Tune Plus window, observe the mass spectra of the singly-charged ions of calibration solution. The ions are as follows. See Figure 22.
 - Caffeine: *m/z* 195
 - MRFA: *m/z* 524
 - Ultramark 1621: *m/z* 1022, 1122, 1222, 1322, 1422, 1522, 1622, 1722, 1822

Note Based on the LC flow rate of your experiment, you can specify the value of each of the following tuning parameters on the LTQ XL MS detector: sheath, Auxiliary, and Sweep gas pressures, ESI needle (or "spray") voltage, ion transfer capillary temperature, and probe position. Automatic tuning sets the values of the other parameters.

- b. At the top of the Spectrum view, notice the values for the ionization time (IT) and normalization level (NL). See Figure 22.
- c. Click the **API Source** button to open the ESI Source dialog box. (See the Spray Current readback shown in Figure 19.)



Testing the Operation of the MS Detector in the ESI/MS Mode

- d. Observe the value for the Spray Current readback and the values for *NL* and *IT* in the Spectrum view. As calibration solution infuses, and the readback values fluctuate, ask yourself the following questions about the ion current signal:
- Is the signal present?
- Is the signal stable, varying by less than about 15% from scan to scan?

If you answered "yes" to the questions in step 4.d, then your MS detector is operating properly.

If you answered "no" to either of these questions, try the following troubleshooting measures:

- Ensure that the fused-silica sample tube does not extend beyond the tip of the ESI needle.
- Ensure that the entrance to the ion transfer capillary is clean, and is not covered with a piece of septum.
- Ensure that the solution entering the probe is free of air bubbles and that the tubing and connectors are free of leaks.

Congratulations! You have demonstrated that your MS detector is operating properly in the ESI mode. You are now ready to tune and calibrate the MS detector. Leave your LTQ XL MS detector as it is, and go to the next section, Tuning the MS Detector Automatically in the ESI/MS Mode.



Figure 22. Spectrum view of the Tune Plus window, showing ionization time (IT) and normalization level (NL) of the calibration solution

Tuning the MS Detector Automatically in the ESI/MS Mode



You tune the MS detector automatically in the ESI/MS mode to optimize important parameters, including heated capillary voltage and tube lens voltage.

Tune the MS detector automatically

- 1. On the Control/Scan Mode toolbar, click the **Tune** button to display the Tune dialog box.
- 2. If necessary, click the **Automatic** tab to display the Automatic tuning page. See Figure 23.
- 3. In the What to Optimize On group box, select the Mass option button to make active the Mass spin box.

Tune	×
Automatic Semi-Automatic Manual Collision Energy	
What to Optimize On	
O Base Peak	
• Mass (m/z): 195.10	
- Status	
Start Cancel Print Help	

Figure 23. Tune dialog box, showing the Automatic tuning page

4. In the Mass spin box, enter **195.1** to specify that the LTQ XL MS detector optimize your Tune Method on the peak at *m*/*z* 195.1.

Note In this example, you use the mass peak at m/z 195.1 to optimize the Tune Method. However, you can optimize the tune on any mass peak of the calibration solution.

- 5. Start the automatic tuning procedure, as follows:
 - a. Click **Start**. A message box displays the following message:

Please ensure that the 500 microliter syringe is full.

Ensure that the syringe contains at least 450 μ L calibration solution.

- b. Click **OK** to close the message box, and return to the Tune dialog box.
- 6. On the File/Display toolbar, click the **Graph View** button to display the Graph view. See Figure 24.
- 7. Observe the Tune Plus window and the Tune dialog box. While automatic tuning is in progress, the LTQ XL MS detector displays various tests in the Spectrum and Graph views in Tune Plus and displays various messages in the Status group box in the Tune dialog box. Your Tune Plus window should now look similar to the one shown in Figure 24.
- 8. Click the **ESI Source** dialog box to examine the ESI source parameters after tuning. Compare the settings shown in Figure 25 with the pre-tune settings shown in Figure 19 on page 3-49.
- 9. Click the **Ion Optics** toolbar button to display the Ion Optics dialog box. The parameters in the Ion Optics dialog box are optimized automatically by the LTQ XL MS detector. See Figure 26.

You have now successfully tuned the MS detector in ESI/MS mode using the calibration solution. Go to the next section, Saving Your ESI/MS Tune Method.



3 Tuning and Calibrating Automatically in the ESI/MS Mode

Tuning the MS Detector Automatically in the ESI/MS Mode



Figure 24. Tune Plus window, showing the results of a typical automatic tune procedure

ESI Source	×
	Actual
Sheath Gas Flow Rate (arb): 10	9.90
Aux <u>G</u> as Flow Rate (arb): 0	0.01
S <u>w</u> eep Gas Flow Rate (arb): 0	0.00
Spray Voltage (kV) : 4.50	4.50
Spray Current (μΑ):	0.05
Capillary Temp (*C): 275.00	274.99
Capillary Voltage (V): 27.00 📫	27.04
<u>I</u> ube Lens Offset (V): 100.00 📑	100.15
Apply OK Cancel <u>H</u> e	elp



Ion Optics	×
	Actual
Multipole 00 Offset (V): -5.50	-5.51
Intermultipole Lens 0 Voltage (V): -5.50	-5.50
Multipole 0 Offset (V): -5.25	-5.24
Intermultipole Lens 1 Voltage (V): -10.00	-9.98
Gate Lens Voltage (V): -52.00	-52.00
Multipole 1 Offset (V): -6.50	-6.50
Multipole <u>B</u> F Amplitude (V p-p): 500.00	500.02
Eront Lens (V): -6.50	-6.50
Apply OK Cancel H	łelp
Apply OK Cancel E	

Figure 26. Ion Optics dialog box, showing examples of voltages of lenses and Intermultipoles, which are optimized by the LTQ XL automatic tuning procedure

Saving Your ESI/MS Tune Method

You can save the parameters you just set in a Tune Method specific to your particular analyte and solvent flow rate. (In this case, you save settings obtained using calibration solution.) You can recall the Tune Method and use it as a starting point for optimizing the MS detector on a different analyte of interest or at a different flow rate.

Note You must save the Tune Method while the MS detector is On.

Save your ESI/MS Tune Method (for low-flow operation) when automatic tuning is complete

1. Choose File > Save As to display the Save As dialog box. See Figure 27.

Save As			<u>?</u> ×
Save in: 🗀	methods	▼ ← 🗈	💣 🎟 •
APCI.LTQT APCIHighFl AutoTune.L APCILowFl APCITune.I AutoTune.L CALI.LTQT	une ow.LTQTune .TQTune ow.LTQTune LTQTune .TQTune une	 ESI.LTQTune ESIHighFlow.LTQTune ESILowFlow.LTQTune ESITune.LTQTune 	
File <u>n</u> ame:	ESImyTune		<u>S</u> ave
Save as <u>t</u> ype:	Tune Files (*.LTC]Tune)	Cancel
No file select	mation		



2. Select the C:\Xcalibur\methods folder.
- 3. Click the **File Name** text box, and type **ESImyTune** to name the Tune Method ESImyTune.LTQTune.
- 4. Click **Save** to save the Tune Method, and return to the Tune Plus window. Note that the Tune Method is named ESImyTune.LTQTune.

Once you have tuned the MS detector, you are now ready to calibrate.

Calibrating the MS Detector Automatically

Calibrate the MS detector automatically from the Tune Plus window

- 1. Choose **Control > Calibrate** to display the Calibrate dialog box.
- 2. If necessary, click the **Automatic** tab to display the Automatic calibration page. See Figure 28.

Calibrate	X
Automatic Semi-Automatic Check	
Calibration Items	
Multipole RF Frequency	
Main RF Frequency	
Electron Multiplier Gain	
Mass and Resolution for Normal Scan Types	
Mass and Resolution for ZoomScan and UltraZoom Types	
Isolation and Activation Waveforms	
Status	
· · · · · · · · · · · · · · · · · · ·	
Set Instrument to Standby when <u>Finished</u>	
Start Cancel Print Help	



- 3. Start the automatic calibration procedure, as follows:
 - a. Click Start. A message box displays the following message:
 Please ensure that the 500 microliter syringe is full.

Ensure the syringe contains at least 450 µL calibration solution.

- b. Click **OK** to close the message box, and return to the Calibrate dialog box.
- 4. Observe the Tune Plus window and the Calibrate dialog box. While the automatic calibration is in progress, the LTQ XL MS detector displays a variety of test results in the Spectrum and Graph views and displays a variety of messages in the Status box of the Calibrate dialog box.

The automatic calibration procedure typically takes about 40 min.

When the LTQ XL MS detector completes the calibration procedure it restores the full scan ESI mass spectrum in the Spectrum view. The Instrument Messages dialog box is displayed, which indicates whether or not the calibration procedure for an item is successful.

- If a calibration item is successful, the LTQ XL MS detector saves the new calibration parameter automatically to the hard disk.
- If a calibration item fails, you can try calibrating on that item again after you ensure the following: the spray is stable, the solution flow rate is sufficient, and all the ions in the calibration solution are present with adequate signal-to-noise ratios. If the sensitivity of the ions is low, increase the solution flow rate somewhat, and then use the *semi-automatic* calibration procedure to calibrate the specific parameter that failed. See Figure 29. Consider deselecting the ZoomScan Mode option if repeated failures occur.

When all calibration items are successful, your MS detector is properly tuned and calibrated for low-flow experiments. A successful calibration exhibits adequate intensities of the following calibrant ions: m/z 195, 524, 1222, 1522, and 1822. In many cases, fine tuning on your particular analyte is not necessary if the intensity of these ions is sufficient. You are ready to analyze samples if you do not need to maximize the intensity of the ion signals for your analyte.

The procedures for changing the solution flow rate and optimization of the MS detector parameters for reserpine, or your particular analyte, are explained in Chapter 4: "Tuning with Your Analyte in LC/ESI/MS Mode". Before you tune with your analyte, go to the next section, Cleaning the MS Detector after Tuning and Calibrating.

3 Tuning and Calibrating Automatically in the ESI/MS Mode

Calibrating the MS Detector Automatically



Figure 29. Tune Plus window with Calibrate dialog box, showing the results of a successful semi-automatic calibration procedure

3 Tuning and Calibrating Automatically in the ESI/MS Mode Cleaning the MS Detector after Tuning and Calibrating

Cleaning the MS Detector after Tuning and Calibrating



This topic describes how to clean your MS detector after using the calibration solution, in preparation for acquiring data on your analyte of interest.

Clean the MS detector after calibrating

1. Click **On/Standby** to put the MS detector in Standby mode. When the MS detector is in Standby, the LTQ XL MS detector turns off the sheath gas, Auxiliary gas, Sweep gas, ESI high voltage, and syringe pump. The MS detector stops scanning, and the LTQ XL MS detector freezes the displays for the Spectrum and Graph views.



CAUTION Always place the MS detector in Standby (or Off) before you open the API source to atmospheric oxygen. The presence of oxygen in the ion source when the MS detector is On could be unsafe. (The LTQ XL MS detector automatically turns the MS detector Off when you open the API source, however, it is best to take this added precaution.

- 2. Remove the syringe from the syringe pump holder, as follows:
 - a. Squeeze the blue buttons, and pull back on the syringe pump handle to free the syringe.
 - b. Remove the syringe from the holder.
 - c. Disconnect the tip of the syringe needle from the Teflon tubing.
- 3. Clean the syringe thoroughly, as follows:
 - a. Clean the syringe with a solution of 5% formic acid in water.
 - b. Rinse the syringe with a solution of 50:50 methanol:water.
 - c. Use acetone to rinse the syringe. (Repeat this step several times.)
- 4. To gain access to the ion transfer capillary, the Ion Max ion source housing and the ion sweep cone need to be removed. Refer to the topic "Removing the Ion Max Ion Source Housing" on page 33 for instructions for removing the Ion Max ion source housing.
- 5. Remove the ion sweep cone as follows:
 - a. Put on a pair of talc-free gloves.

3 Tuning and Calibrating Automatically in the ESI/MS Mode Cleaning the MS Detector after Tuning and Calibrating



CAUTION AVOID BURNS. At operating temperatures, the ion transfer tube can severely burn you! The ion transfer tube typically operates between 200 and 400 °C. Always allow the ion sweep cone to cool to room temperature (for approximately 20 min) before you touch or remove this component. Always be careful not to touch the entrance end of the ion transfer tube when it is exposed.

- b. Grasp the outer ridges of the ion sweep cone and pull the cone straight off of the API cone seal. Note, you might need to loosen the set screws on the ion sweep cone in order to remove it.
- 6. Remove the ion transfer capillary by using the custom tool provided.
- 7. Clean the ion sweep cone and the ion transfer capillary as follows:
 - a. Place the ion sweep cone and the ion capillary tube in a beaker of 50:50 methanol/water.
 - b. Sonicate the components for 15 min.
- 8. Reinstall the ion transfer capillary.
- 9. Reinstall the ion sweep cone as described in "Installing the Ion Sweep Cone" on page 34.
- 10. Place a small Teflon-coated septum over the entrance end of the ion transfer capillary to seal the vacuum chamber of the MS detector.
- 11. Flush the sample transfer line, sample tube, and ESI probe thoroughly with a solution of 5% formic acid in water (or with another appropriate solvent), as follows:

Note The solvent that you use to flush the sample transfer line, sample tube, and ESI probe assembly depends on the solvent system you use to dissolve your samples. For example, if you are using a buffered solution of a high concentration, an acidic solution is appropriate.

- a. Fill a clean, 250 μ L Unimetrics syringe with a solution an appropriate solvent.
- b. While holding the plunger of the syringe in place, carefully insert the needle of the syringe into the free end of the Teflon tube.
- c. Flush the sample transfer line, sample tube, and ESI probe with the solution by slowly depressing the syringe plunger. Visually confirm that the solution is exiting the tip of the ESI probe on the inside of

the probe assembly. Use a lint-free tissue to gently remove the excess solution as it exits the probe.

- d. Remove the needle of the syringe from the Teflon tube.
- 12. Repeat step 11 with a solution of 50:50 methanol:water.
- 13. Repeat step 11 with acetone.
- 14. Clean the spray shield as follows:
 - a. Fill a spray bottle with solvent solution.
 - b. Temporarily place a large Kimwipe (or other lint-free tissue) beneath of the spray shield. (The Kimwipe is required to absorb the solution used to flush the ion transfer capillary and spray shield.)
 - c. Use the spray bottle to flush contaminants from the exterior surface of the spray shield.
 - d. Remove the Kimwipe you used to absorb the solution. Swab the surface of the spray shield with a dry Kimwipe.
 - e. Repeat step 14.a through step 14.d with acetone to remove the (high molecular weight) Ultramark 1621.
- 15. Being careful not to touch the ion transfer capillary with your hand, remove the septum from the entrance end of the ion transfer capillary.
- 16. Reinstall the Ion Max ion source housing as described in "Installing the Ion Max Ion Source Housing" on page 36.

The MS detector is now clean and ready for acquiring data on your analyte of interest.

If you plan to run analytical samples in high-flow ESI mode (using flow rates between 50 and 1000 μ L/min), the procedures in Chapter 4: "Tuning with Your Analyte in LC/ESI/MS Mode" explain how to optimize the tune for this situation.

Chapter 4 Tuning with Your Analyte in LC/ESI/MS Mode

This chapter provides information on tuning the MS detector in the LC/ESI/MS mode using your analyte. You optimize the sensitivity of the MS detector to your analyte through an automatic procedure.

The customized Tune Methods contained in your LTQ XL data system are optimized for a wide range of applications, and they can be used without further tuning of your MS detector. However, for certain applications you might need to tune and optimize several MS detector parameters.

For instance, the most important parameters that interact with the ESI interface and signal quality are as follows:

- Electrospray voltage
- Heated capillary temperature (voltage)
- Tube lens voltage
- Capillary voltage
- Sheath gas flow rate
- Auxiliary gas flow rate
- Sweep gas flow rate

The settings for these parameters depend on the solvent flow rate and target analyte composition. In general, you should fine tune your MS detector whenever you change the solvent flow rate conditions of your particular application. In this procedure, you use the ESI low-flow Tune Method *ESImyTune.LTQTune* as a starting point, then further optimize the MS detector parameters using an automatic procedure. The automatic procedure adjusts the tube lens voltage, capillary voltage, and voltages applied to the ion optics until the ion transmission of your analyte is maximized.

The capillary is heated to maximize the ion transmission to the MS detector. For ESI only, you set the ion transfer capillary temperature proportional to the flow rate of your solution. Refer to Table 1-2 for guidelines for setting operating parameters for LC/ESI/MS. For this procedure, the ion transfer capillary temperature is set to 350 °C, and the sheath gas is set to 30.

Note

- 1. If your experiment is performed at a flow rate below 10 μ L/min, and the results you want can be obtained without optimizing the MS detector on your particular analyte, go to Chapter 5: "Acquiring ESI Sample Data Using the Tune Plus Window" to acquire sample data.
- 2. Before you optimize the tune for your analyte of interest, ensure that the LTQ XL MS detector has been calibrated within the previous three months. If the system needs to be calibrated, refer to the procedures in Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode".

To tune the MS detector in the ESI/MS (high-flow) mode using your analyte, perform the following tasks:

- Set up the MS detector for your specific analyte from the Tune Plus window.
- Infuse your analyte into the MS detector using a syringe pump connected to the LC with a Tee union.
- Optimize the MS detector parameters for your analyte while the solution flows into the MS detector.

This chapter contains the following topics:

- Setting Up to Introduce Sample by Syringe Pump into Solvent Flow from an LC
- Setting Up to Tune the MS Detector with Your Analyte
- Optimizing the MS Detector Tune Automatically with Your Analyte
- Saving the ESI/MS Tune Method

Setting Up to Introduce Sample by Syringe Pump into Solvent Flow from an LC

This topic describes setting up the MS detector to introduce your analyte by syringe pump into solvent flow from an LC.

Plumbing connections for ESI/MS sample introduction into LC solvent flow from the syringe pump

- 1. Connect a 4 cm (1.5 in.) segment of Teflon tubing with a (brown) fingertight fitting and a (brown) ferrule to the (black) LC union. (See Figure 17 on page 46.)
- 2. Fill a clean, 500-µL Unimetrics syringe with the 125 fg/µL solution of reserpine or your analyte of interest. (See Appendix A: "Sample Formulations" for a procedure for making the reserpine tuning solution.)
- 3. Insert the needle of the syringe into the segment of Teflon tube. Check that the needle tip of the syringe fits readily into the opening in the free end of the Teflon tubing. If necessary, you can enlarge the opening in the end of the tubing slightly.
- 4. Place the syringe into the syringe holder of the syringe pump.
- 5. While squeezing the blue release buttons on the syringe pump handle, push the handle forward until it just contacts the syringe plunger.
- 6. Connect the fused-silica infusion line from the (black) LC union to the (black) LC Tee union, as follows. See Figure 30.
 - a. Connect the infusion line with a (brown) fingertight fitting and a (brown) ferrule to the free end of the LC union.
 - b. Connect the other end of the infusion line with a (red) fingertight fitting and a (brown) ferrule to the side arm of the LC Tee union.
- 7. Connect an appropriate length of (red) PEEK tubing from the (stainless steel) grounding union to the (black) LC Tee union, as follows:
 - a. Use a PEEK tubing cutter to cut a 4 cm (1.5 in.) length of the PEEK tubing.
 - b. Connect the PEEK tubing with a (brown) fingertight fitting and a (brown) ferrule to the grounding union.
 - c. Connect the other end of the PEEK tubing with a (brown) fingertight fitting and a (brown) ferrule to the LC Tee union.

Setting Up to Introduce Sample by Syringe Pump into Solvent Flow from an LC



Figure 30. ESI/MS plumbing connections for the LC Tee union

8. If you have not already done so, connect the PEEK safety sleeve and fused-silica sample tube from the grounding union to the ESI probe sample inlet as described in the section, *Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve* in the *Ion Max API Source Hardware Manual*.

If you have installed the stainless steel needle in the ESI probe, connect the PEEK safety sleeve and fused-silica capillary tube from the grounding union to the ESI probe sample inlet as described in the section, *Installing a New Stainless Steel Needle in the ESI Probe* and *Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve* in the *Ion Max API Source Hardware Manual.*

- 9. Connect an appropriate length of PEEK tubing (transfer line from the divert/inject valve) from the divert/inject valve to the free end of the (black) LC Tee union, as follows.
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and a (stainless steel) ferrule to port 3 of the divert/inject valve. See Figure 31.

b. Connect the other end of the PEEK tubing with a (brown) fingertight fitting and a (brown) ferrule to the free end of the LC Tee union. (See Figure 30.)



- **Figure 31.** Divert/Inject valve, showing the correct setup for tuning by syringe infusion and showing the flow of liquid through the valve in the Detector and Waste positions
- 10. Connect an appropriate length of PEEK tubing (transfer line from the LC) from the divert/inject valve to the LC, as follows:
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and a (stainless steel) ferrule to port 2 of the divert/inject valve.
 - b. Connect the other end of the PEEK tubing with a proper fitting and a ferrule to the outlet of the LC.
- 11. Connect an appropriate length of PEEK tubing (waste line) from the divert/inject valve to a waste container, as follows:
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and a (stainless steel) ferrule to port 1 of the divert/inject valve.
 - b. Insert the other end of the PEEK tubing in a suitable waste container.

The MS detector is now properly set up to introduce your analyte by syringe pump into solvent flow from an LC.

Setting Up to Tune the MS Detector with Your Analyte



Set up the MS detector to tune automatically on your analyte in ESI/MS mode

In this procedure, you can use the reserpine solution described in Appendix A: "Sample Formulations", or you can use a solution of an analyte of interest to you.

CAUTION Do not infuse calibration solution at flow rates above 10 μ L/min. Ultramark 1621 can contaminate your system at high concentrations.

Note The following procedures assume that you are familiar with your LTQ XL instrument and the Tune Plus window. If you need additional guidance, see LTQ XL online Help and/or the *LTQ XL Hardware Manual*.

- 1. Open the Tune Plus window from the Start button on your Windows XP Desktop, as follows:
 - a. Choose **Start > Programs > Xcalibur > Xcalibur** to display the Xcalibur Home Page Roadmap view.
 - b. Click the Instrument Setup button to display the Instrument Setup window.
 - c. Click the LTQ XL button to display the New Method page.
 - d. Click the Tune Plus button to display the Tune Plus window.
- 2. In Tune Plus, click the **On/Standby** button to take the MS detector out of Standby mode and turn it On. The MS detector begins scanning, the LTQ XL MS detector applies high voltage to the ESI probe, and the LTQ XL MS detector shows a real-time display in the Spectrum view.
- 3. Open the ESImyTune.LTQTune Tune Method, the Tune Method you saved in Chapter 3, as follows:
 - a. On the File/Display toolbar, click the **Open File** icon to display the Open dialog box.
 - b. Browse to the folder C:\Xcalibur\methods. Then, select the file ESImyTune.LTQTune.
 - c. Click **Open** to open the file. Tune Plus downloads the Tune Method parameters to the MS detector, and the title bar in the Tune Plus window should read as follows:
 C:\Xcalibur\methods\ESImyTune.LTQTune Tune Plus





4. Define the scan parameters for tuning with your analyte in the ESI/MS mode, as follows:



- a. In the Instrument Control toolbar, click **Define Scan** to open the Define Scan dialog box. See Figure 32.
- b. In the Scan Description group box, in the Mass Range list box, select *Normal* to allow for a selection of mass ranges between m/z 150 to 2000.
- c. In the Scan Rate list box, select *Normal* to specify a normal scan rate.
- d. In the Scan Type list box, select *SIM* specify a Selected Ion Monitoring experiment.
- e. In the Scan Time group box, in the Number of Microscans spin box, type 1 to set the total number of microscans to 1.
- f. In the Max. Inject Time spin box, type *200.000* to specify a 200 ms maximum injection time.
- g. In the Scan Ranges group box, in the Input list box, select Center/Width to make available the Center Mass and Width text boxes in the Scan Ranges table.
- h. In the Source Fragmentation group box, confirm that the On check box is not selected (
) to specify that the ion source fragmentation option is turned off.
- i. In the Scan Ranges group box (Scan Ranges table), in the Center Mass text box, type the mass of your analyte to set the center of mass for the scan range. If reserpine is the analyte you would enter **609.20** to set the center mass for the scan range to m/z 609.20. If you are using another analyte enter its mass value into the Center Mass text box (Figure 32)
- j. In the Width text box, type **2.0** to set the width of the scan range to 2.0 daltons.
- k. Verify that the settings in your Define Scan dialog box are the same as those shown in Figure 32.
- 1. Click **OK** to apply the MS detector scan parameters and to close the Define Scan dialog box.

Setting Up to Tune the MS Detector with Your Analyte

Define Scan									×
Scan <u>H</u> istory: ITMS + c SIM ms [60	08.20-6	10.20]							- 🖻 🛍
Scan Description	_ MSr	n Settings					Sca	an Ranges —	
Mass Range: Normal 💌	n	Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)	#	Center Mass (m/z)	Width (m/z)
	2		1.0	20.0	0.250	10.000	1	609.20	2.00
Scan Lype: SIM	3		1.0	20.0	0.250	10.000	2		
	4		1.0	20.0	0.250	10.000	3		
- Soon Time	5		1.0	20.0	0.250	10.000	4		
	- 5		1.0	20.0	0.250	10.000	<u> </u>		
	<u>-</u>		1.0	20.0	0.250	10.000			
Max. Inject Time (ms): 200.000			1.0	20.0	0.250	10.000			
	10		1.0	20.0	0.250	10.000			
Source Fragmentation			tivation	2010	0.200	,0.00	10 10	nput: Center	Width 💌
Apply OK Cancel Help Injection RE Activation									

Figure 32. Define Scan dialog box, showing typical settings for acquiring reserpine data of the SIM type

١Ļ

- 5. On the Control/Scan Mode toolbar, click **Centroid/Profile** to toggle the data type to profile. (The picture on the button should be the same as that shown here.)
- \odot
- 6. Click **Positive/Negative** to toggle the ion polarity mode to positive. (The picture on the button should be the same as that shown here).

You have completed setting up to tune your MS detector with your analyte in ESI/MS mode.

Optimizing the MS Detector Tune Automatically with Your Analyte

Ų



Optimize the MS detector tune automatically to maximize the ion transmission of reserpine (or your analyte of interest) for a high-flow experiment. Thermo Electron recommends that you begin optimizing after you have successfully passed an automatic tuning procedure and an automatic calibration procedure with the calibration solution infused at 5 μ L/min.

The following procedure describes how to optimize the MS detector Tune Method with the reserpine m/z 609.2 peak at an LC flow rate of 400 μ L/min. You can also carry out this procedure with your analyte of interest and at your particular LC flow rate. (Refer to Table 2 for guidelines about setting flow rates and temperatures.)

Optimize the MS detector Tune Method

- 1. On the Control/Scan Mode toolbar, click the **Tune** button to display the Tune dialog box.
- 2. If necessary, click the **Automatic** tab to display the Automatic tuning page. See Figure 33.
- 3. In the What to Optimize On group box, select the Mass option button to make active the Mass spin box.
- 4. In the Mass spin box, enter **609.2** (or the appropriate mass of your analyte of interest) to specify that the LTQ XL MS detector is to use the peak at m/z 609.2 (or the appropriate mass of your analyte of interest) to optimize your tune.
- 5. Ensure that the Divert/Inject valve is in the Detector position, as follows:
 - a. Click the **Divert/Inject** button to open the Divert/Inject Valve dialog box. See Figure 34.
 - b. Click the **Detector** option button.
 - c. Click Close.
- 6. Start the automatic tuning procedure from the Tune dialog box, as follows:
 - a. Click **Start**. A message box displays the following message:

Please ensure that the 500 microliter syringe is full.

Ensure the syringe pump contains at least 450 μL of the 125 fg/ μL reserpine tuning solution.

- b. Click **OK** to close the message box, and return to the Tune Plus window.
- 7. In the File/Display toolbar, click the **Graph View** button to display the Graph view.

Tune X
Automatic Semi-Automatic Manual Collision Energy What to Optimize On Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Collision Energy Image: Collision Energy Image: Image: Energy Image: Energy Image: Collision Energy Image: Image: Energy Image: Energy Image: Collision Energy Image: Image: Energy Image: Energy Image: Energy Image: Image: Energy Image: Energy Image: Energy Image: Image: Energy Image: Energy Image: Energy Image: Image: Image: Image: Energy Image: Energy Image: Energy
Status
<u>S</u> tart Cancel <u>Print</u> <u>H</u> elp

Figure 33. Tune dialog box, showing the Automatic tuning page



Divert/Inject V	×	
Load	Inject	
۲	0	
Detector	Waste	
Close	<u>H</u> elp	
		-

Figure 34. Divert/Inject Valve dialog box

8. Observe the Tune Plus window and the Tune dialog box. While automatic tuning is in progress, the LTQ XL MS detector displays various tests in the Spectrum and Graph views in the Tune Plus window and displays various messages in the Status group box in the Tune dialog box. Your Tune Plus window should now look similar to the one shown in Figure 35.

Note The most important parameters that affect the signal quality during ESI/MS operation are electrospray voltage, ion transfer capillary temperature, heated capillary voltage, tube lens voltage, gases, and solution flow rate. If any one of these parameters is changed, you need to reoptimize MS detector parameters. You can use the Semi-Automatic tune procedure to tune the MS detector on individual parameters.

You have now successfully tuned the MS detector in ESI/MS mode for the compound reserpine (or your analyte of interest).

4 Tuning with Your Analyte in LC/ESI/MS Mode

Optimizing the MS Detector Tune Automatically with Your Analyte



Figure 35. Tune Plus window with the Tune dialog box, showing the Automatic tuning page

Saving the ESI/MS Tune Method

Save your ESI/MS Tune Method (for a high-flow experiment using your analyte) when automatic tuning is complete

Note Save the Tune Method while the MS detector is On, if any of the ion source parameters are different from those with which you started.

1. Choose **File > Save As** to display the Save As dialog box. See Figure 36.

Save As			? ×
Save jn: 🗀	methods	- + 1	📸 🎟 -
🥼 Alprazolam.	LTQTune		
APCITune.	LTQTune TQTune		
ESITune.L	TQTune		
File <u>n</u> ame:	reserpine		<u>S</u> ave
Save as <u>t</u> ype:	Tune Files (*.LTQTune)	•	Cancel
- Header Infor	mation		
No file selec	sted		
			//



- 2. Select the C:\Xcalibur\methods folder.
- 3. Click the **File Name** text box, and enter **reserpine** (or the name of your analyte of interest).
- 4. Click **Save** to save the Tune Method, and return to the Tune Plus window. Note that the Tune Method is named reserpine.LTQTune.

The Tune Method is now properly saved and you are ready to acquire data on your analyte of interest.

Chapter 5 Acquiring ESI Sample Data Using the Tune Plus Window

This chapter describes how to acquire LC/ESI/MS sample data using the Tune Plus window. This experiment uses reserpine but you can use the same procedure with your analyte of interest.

Note The following procedures assume that you are familiar with your LTQ XL instrument and the Tune Plus window. If you need information, refer to the LTQ XL online Help, *LTQ XL Getting Connected*, and/or the *LTQ XL Hardware Manual*.

Ensure that you have completed the procedures in Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode" and Chapter 4: "Tuning with Your Analyte in LC/ESI/MS Mode".

This chapter contains the following sections:

- Setting Up to Acquire MS/MS Data in the Full Scan Type
- Setting Up to Introduce Sample by Loop Injection into Solvent Flow from an LC
- Acquiring MS Data in the SIM Scan

Setting Up to Acquire MS/MS Data in the Full Scan Type

Prepare to acquire MS/MS data in the Full scan type on reserpine (or on your analyte of interest). You need to optimize the isolation width and the relative collision energy parameters before you acquire MS/MS data.

You first optimize the isolation width to ensure that the ion of interest is isolated effectively, and then you optimize the collision energy to ensure that fragmentation of the parent ion is efficient. The relative collision energy for a particular analysis depends on the type of sample you are analyzing.

The information in this topic applies to operation of the LTQ XL MS detector in either the ESI or the APCI mode.

This topic contains the following subtopics:

- Optimizing the Isolation Width and Setting Up to Optimize the Collision Energy
- Optimizing the Collision Energy Automatically for an MS/MS Experiment

Optimizing the Isolation Width and Setting Up to Optimize the Collision Energy





- data for rese
- ⇒ ●

Optimize the isolation width and set up to optimize the collision energy for an MS/MS experiment

Note The collision energy is optimized on the LTQ XL MS detector by changing the values for the parameter Normalized Collision Energy in the MSⁿ Settings group box of the Define Scan dialog box. For this experiment, and for most applications, leave the parameters Activation Q and Activation Time set to their default values. For more information about these parameters, refer to the online Help.

- 1. If you have not already done so, from the Tune Plus window, click the **On/Standby** button to take the MS detector out of Standby mode and turn it On.
- 2. Ensure that the Centroid data type is selected. (The picture on the button should be the same as that shown here.)
- 3. Ensure that the scan parameters are defined to acquire MS/MS Full scan data for reserpine (or your analyte of interest), as follows:
 - a. Click the Define Scan button to open the Define Scan dialog box. See Figure 37.

Define Scan											×
Scan <u>H</u> istory: ITMS + c Full ms [16	5.00-6	50.00]								- 🖻 🛍	
Scan Description	_ MSr	n Settings						Sca	n Ranges—		1
Mass Range: Normal 💌	n	Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)		#	First Mass (m/z)	Last Mass (m/z)	
	2	609.20	3.0	20.0	0.250	10.000		1	165.00	650.00	
Scan <u>Type</u> : Full	3		1.0	20.0	0.250	10.000		2			
	4		1.0	20.0	0.250	10.000		3			
	5		1.0	20.0	0.250	10.000		4			
Scan Time	6		1.0	20.0	0.250	10.000		5			
Mi <u>c</u> roscans: 1	7		1.0	20.0	0.250	10.000		6			
Max Inject Time (ms): 200.000	8		1.0	20.0	0.250	10.000		- 7			
	9		1.0	20.0	0.250	10.000		8			
	10		1.0	20.0	0.250	10.000		9			
Source Fragmentation								10			
□ <u>0</u> n <u>Energy</u> (V): 20.0 <u>+</u>		<u>W</u> ideband Ac	tivation					Ī	nput: From/	To 💌	
	Арр	y _	OK .	Cancel	Hel	p	In	jectic	n R <u>E</u>	Acti <u>v</u> ation	

- **Figure 37.** Define Scan dialog box, showing initial settings to optimize the isolation width of an MS/MS experiment for reserpine
 - b. Verify that the values in your dialog box are the same as those shown in Figure 37. Start with a relatively wide Isolation Width. Leave the Define Scan dialog box open.
 - 4. At this time you might want to turn on your LC pump and specify a flow rate of 0.4 mL/min, for example, to ensure that your system does not leak.



- 5. Click the **Syringe Pump** button to display the Syringe Pump dialog box. See Figure 38.
- 6. Turn on the syringe pump and set an infusion flow rate of 5 $\mu L/\text{min},$ as follows:
 - a. In the Flow Control group box, click the **On** option button to make active the Flow Rate spin box.
 - b. Type 5 in the Flow Rate spin box to specify a rate of 5 μ L/min.

Note This procedure assumes that you are using the $500-\mu L$ Unimetrics syringe that is provided with your LTQ XL system. If you are using another type of syringe, select the option button corresponding to your syringe. Setting Up to Acquire MS/MS Data in the Full Scan Type

S	vringe Pumn				>	٢l
	- Flow Control					-
					Actual	
	⊙ <u>O</u> n	Flow <u>R</u> ate (μ L/min):	5.00	·	0.00	
	O Ołł					
	Туре]	
	⊖ Ha <u>m</u> ilton	⊻olume (μL):	500	•		
	• <u>U</u> nimetrics	Syringe [D (mm):	3.260	* *		
	O Oth <u>e</u> r					
	Apply	ОК	Cancel	<u>⊦</u>	lelp	

Figure 38. Syringe Pump dialog box

- c. If you are using a standard Unimetrics or Hamilton syringe, set up the syringe parameters as follows:
 - i. In the Type group box, click the **Unimetrics** or **Hamilton** option button to specify the proper syringe type.
 - ii. Click the **Volume** list box arrow to display the list of available volumes, and then select *500* (or your syringe size) from the list to set the proper syringe volume. Note that, if you are using a Unimetrics syringe, the LTQ XL MS detector automatically sets the syringe ID to its proper value of 3.260 mm.
- d. If you are not using a Unimetrics or Hamilton syringe, set up the syringe parameters as follows:
 - i. In the Type group box, click the **Other** option button to make active the syringe ID spin box.
 - ii. Enter the inner diameter of your syringe in the Syringe ID spin box.
- e. Click **Apply** to apply the syringe parameters and start the syringe pump.
- f. Finally, move the Syringe Pump dialog box out of the way, to the top of the monitor screen.
- 7. In the Tune Plus window, observe the mass spectrum of reserpine (or your analyte of interest). Also observe the values for NL and IT

(*Normalization Level* and *Ion Time*), while you optimize the value of the Isolation Width in the Define Scan dialog box, as follows:

- a. In the Define Scan dialog box, in the MS^n Settings group box, in the Isolation Width box, type 3 to specify an isolation width of m/z 3. Then, click **Apply**.
- b. In the Tune Plus window, observe the mass spectrum for the parent ion of reserpine, m/z 609.2. Ensure that the readback values for NL and IT are relatively stable.
- c. Again, in the Define Scan dialog box, in the MS^n Settings group box, in the Isolation Width box, type **2.8** to specify an isolation width of m/z 2.8. Then, click **Apply**.

Note The optimum value for the Isolation Width is the smallest m/z width (instrument minimum width =m/z 0.4) that gives a mass spectrum of maximum intensity for only the ions of interest. When the optimum Isolation Width is obtained the values for NL and IT are stable and the mass peak for the parent ion is at its maximum intensity and appears symmetrical. An Isolation Width value that is less than the optimum value causes a substaitial drop in the NL reading. A significant drop in sensitivity indicates that the ions of interest are not effectively isolated.

d. Repeat steps b and c above, entering successively smaller values for Isolation Width. Continue to observe the intensity of the mass spectrum of the parent ion, and ensure that the values for *NL* and *IT* are stable with each change you make to the Isolation Width.

Note After the Isolation Width is optimized, you can compensate for minor changes in tune stability by increasing the Isolation Width value a small amount. This adjustment should be no larger than m/z=1.

- 8. In the Define Scan dialog box, in the MSⁿ Settings group box, in the Normalized Collision Energy box, type **20** to specify an initial value of 20 for the collision energy. Click **Apply**.
- 9. In the Tune Plus window, observe the mass spectrum of the product ions of reserpine (or your analyte of interest). If necessary, increase the value for the Normalized Collision Energy in increments of 5% to cause the clear display of product ion mass spectrum. (After each change in value, click **Apply** to implement the change.) See Figure 39.

5 Acquiring ESI Sample Data Using the Tune Plus Window

Setting Up to Acquire MS/MS Data in the Full Scan Type



- 10. When you have clearly identified a product ion mass-to-charge ratio for reserpine (or your analyte of interest), click the **Tune** button to display the Tune dialog box.
- 11. In the Tune dialog box, click the **Collision Energy** tab to display the page. See Figure 40.

Define Scan											×
Scan <u>H</u> istory: ITMS + c Full ms [16	5.00-6	50.00]								- B B	
Scan Description	⊢ MSr	n Settings						Sca	n Ranges—		1
Mass Range: Normal 💌	n	Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)		#	First Mass (m/z)	Last Mass (m/z)	
	2	609.20	2.0	25.0	0.250	10.000		1	165.00	650.00	
Scan <u>I</u> ype: Full 💌	3		1.0	20.0	0.250	10.000		_ 2			
	4		1.0	20.0	0.250	10.000		3			
	5		1.0	20.0	0.250	10.000		4			
Scan Time	6		1.0	20.0	0.250	10.000		5			
Microscans: 1	7		1.0	20.0	0.250	10.000		6			
May Inject Time (me): 200,000	8		1.0	20.0	0.250	10.000		7			
	9		1.0	20.0	0.250	10.000		8			
	10		1.0	20.0	0.250	10.000		9			
Source Fragmentation		□ <u>W</u> ideband Activation						10 _[r	put: From/	To 💌	
Apply OK Cancel Help Injection RE Activation											

Figure 39. Define Scan dialog box, showing typical settings for acquiring MS/MS data in the Full scan type on reserpine

12. Click the **Product Ion Mass** option button to make active the spin box. Type **397.2** to specify the product ion at *m/z* 397.2 for reserpine. The LTQ XL MS detector can optimize collision energy automatically by using this product ion of reserpine.

Tune		×
Automatic Semi-Automatic	Manual Collision Energy	
What to Optimize	What to Optimize On C <u>I</u> IC Product Ion <u>M</u> ass (m/z) <u>397</u>	.20 ÷
	Results Initial Collision Energy: 25 Best Collision Energy: 26	5.00 % 3.00 %
Status		
<u>S</u> tart	Cancel <u>P</u> rint	<u>H</u> elp

Figure 40. Tune dialog box, showing the Collision Energy page

Optimizing the Collision Energy Automatically for an MS/MS Experiment

The optimum relative collision energy is the one that produces the maximum product ion intensity.

Automatically optimize the relative collision energy for the ESI/MS/MS analysis of reserpine (or your analyte of interest)

1. In the Tune dialog box, on the Collision Energy page (Figure 40), click **Start** to start the optimization procedure. A message box displays the following message:

Ensure that the 500 microliter syringe is full.

Ensure the syringe contains at least 450 μL of the 125 fg/ μL reserpine solution.

- 2. Click **OK** to close the message box, and leave the Tune dialog box open. Your Tune Plus window should now look similar to the one shown in Figure 41.
- 3. In the Spectrum view of Tune Plus, observe the MS/MS Full scan spectrum of reserpine (or that of your analyte of interest).
- 4. When the collision energy is optimized, the Accept Optimized Value dialog box appears. See Figure 42.
- 5. Click **Accept** to accept the new collision energy value, and return to Tune Plus. The new value is displayed in the Define Scan dialog box.
- 6. In the Syringe Pump dialog box, select the Off option button to turn off the syringe pump. Click **Close** to close the Syringe Pump dialog box.
- 7. Click **Cancel** to close the Tune dialog box.

After you optimize the relative collision energy, the LTQ XL MS detector is ready to acquire MS/MS data on your analyte of interest.

5 Acquiring ESI Sample Data Using the Tune Plus Window

Setting Up to Acquire MS/MS Data in the Full Scan Type



Figure 41. Tune Plus window, showing MS/MS product ions of reserpine in the Spectrum view



Figure 42. Accept Optimized Value dialog box

Setting Up to Introduce Sample by Loop Injection into Solvent Flow from an LC

Plumbing connections to introduce sample by loop injection into the solvent flow from an LC $\,$

1. If you have not already done so, connect the PEEK safety sleeve and fused-silica sample tube from the grounding union to the ESI probe sample inlet as described in the section *Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve* in the *Ion Max API Source Hardware Manual*.

If you have installed the stainless steel needle in the ESI probe, connect the PEEK safety sleeve and fused-silica capillary tube from the grounding union to the ESI probe sample inlet as described in the section *Installing a New Stainless Steel Needle in the ESI Probe* and *Installing a New Fused-Silica Sample Tube and PEEK Safety Sleeve* in *Ion Max API Source Hardware Manual.*

- 2. Connect an appropriate length of (red) PEEK tubing (transfer line from the divert/inject valve) from the divert/inject valve to the (stainless steel) grounding union, as follows. See Figure 43.
 - a. Connect a length of PEEK tubing fitted with a (stainless steel) nut and a (stainless steel) ferrule to port 3 of the divert/inject valve. See Figure 44.
 - b. Connect the other end of the PEEK tubing with a (brown) fingertight fitting and a (brown) ferrule to the free end of the grounding union.
- 3. Connect a 5 μ L sample loop with (stainless steel) nuts and (stainless steel) ferrules to ports 1 and 4 of the divert/inject valve. (See Figure 44)



Figure 43. ESI/MS plumbing connections for the divert/inject valve and grounding union

Setting Up to Introduce Sample by Loop Injection into Solvent Flow from an LC



- **Figure 44.** Divert/Inject valve, showing the correct setup for analysis by loop injection and showing the flow of liquid through the valve in the Load and Inject positions
- 4. Connect an appropriate length of PEEK tubing (transfer line from the LC) from the divert/inject valve to the LC, as follows:
 - a. Connect a length of the PEEK tubing with a (stainless steel) nut and (stainless steel) ferrule to port 2 of the divert/inject valve.
 - b. Connect the other end of the PEEK tubing with a proper fitting and a ferrule to the outlet of the LC.
- 5. Connect an appropriate length of PEEK tubing (waste line) from the divert/inject valve to a waste container, as follows:
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and (stainless steel) ferrule to port 6 of the divert/inject valve.
 - b. Insert the other end of the PEEK tubing into a suitable waste container.

Acquiring MS Data in the SIM Scan





scan from the Tune Plus window

Acquire a file of reserpine data using the selected ion monitoring (SIM)

Note The LTQ XL MS detector automatically saves the data you acquire on your hard disk.

1. On the Control/Scan Mode toolbar, click **On/Standby** to take the MS detector out of Standby mode and turn it On. The MS detector begins scanning, the LTQ XL MS detector applies high voltage to the ESI probe, and a real-time display shows in the Spectrum view.

- 2. Verify that the Centroid data type is selected. (The picture on the button should be the same as that shown here.)
- 3. Verify that the scan parameters are defined to acquire SIM data for reserpine (or your analyte of interest), as follows:



- a. Click the Define Scan button to open the Define Scan dialog box. See Figure 45.
- b. Compare the values in your dialog box to those in Figure 45 and click **OK**.

Define Scan									ļ
Scan <u>H</u> istory: ITMS + c SIM ms [60	08.20-610.20]								- 🖻 🛍
Scan Description	MSn Settings						-Sca	n Ranges	
Mass Range: Normal 💌	n Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)		#	Center Mass (m/z)	Width (m/z)
	2	1.0	20.0	0.250	10.000		1	609.20	2.00
Scan <u>Type:</u> SIM	3	1.0	20.0	0.250	10.000		2		
	4	1.0	20.0	0.250	10.000		3		
]	5	1.0	20.0	0.250	10.000		4		
Scan Time	6	1.0	20.0	0.250	10.000		5		
Microscans: 1	7	1.0	20.0	0.250	10.000	1	6		
May Initian Time (such 200,000	8	1.0	20.0	0.250	10.000	1	7		
	9	1.0	20.0	0.250	10.000	1	8		
	10	1.0	20.0	0.250	10.000	1	9		
Source Fragmentation	□ <u>W</u> ideband Ac	tivation					10 <u> </u> r	n <mark>put:</mark> Center	Width 💌
Apply OK Cancel Help Injection RE Activation									

Figure 45. Define Scan dialog box, showing typical settings for acquiring reserpine data in the SIM scan type

4. Turn on the LC pump and specify your flow rate of 400 μ L/min. Ensure that your system is free of leaks.

- 5. On the Control/Scan Mode toolbar, click **Acquire Data** to open the Acquire Data dialog box. See Figure 46.
- 6. Specify the acquisition parameters, as follows:
 - a. In the File Name box, enter **reserpine** to specify a filename.
 - b. In the Sample Name box, enter **reserpine** to specify the sample identity. If you are not using reserpine, type the name of your particular analyte.
 - c. Type a comment about your experiment. (For example, describe the scan mode, scan type, ionization mode, sample amount, or method of sample introduction.). The Xcalibur data system includes the comment on hard copies of your data.
 - d. In the Acquire Time box, select the Continuously option button to acquire data continuously (until you stop the acquisition).
- 7. Leave the Acquire Data dialog box open during data acquisition, but move it to a corner of the Tune Plus window.
- 8. Click **Start** in the Acquire Data dialog box to begin acquiring data. The Acquisition Status group box displays the following message.

State: Acquiring

Time (min):



9. Click **Divert/Inject Valve** to open the Divert/Inject Valve dialog box. See Figure 47.

Acquiring MS Data in the SIM Scan

Acquire Data			<u>? ×</u>
Fol <u>d</u> er:	C:\Calibur\Data	Acquire Time	
<u>F</u> ile Name:	reserpine 001	Continuously	<u>S</u> tart
Sample <u>N</u> ame:	reserpine	C Scans 10 -	Pause
<u>C</u> omment:	SIM, APCI, 50pg loop		⊻iew
Instrument <u>M</u> ethod:	Use instrument method C:\Xcalibur\methods\MyMethod.meth	Coto Standby when Finished Acquisition Status State: Idle	Inst. Setup
	Start Mode Immediate O Contact Closure O Divert Valve OK Cancel <u>H</u> elp	Time (min): 0.000	

Figure 46. Acquire Data dialog box, showing typical settings for acquiring data

Divert/Inject V	alve	×
Load	Inject	
0	۲	
Detector	Waste	
Close	<u>H</u> elp	
		-

Figure 47. Divert/Inject Valve dialog box

- 10. Select the Load option button, and overfill the 5- μ L injector loop with the 125 fg/ μ L solution of reserpine (or a solution of your analyte of interest).
- 11. Select the Inject option button to inject the reserpine solution into the ESI source. Leave the Divert/Inject Valve dialog box open.
- 12. Observe the reserpine peak (*m/z* 609.2), or that of your analyte of interest, in the Spectrum view. See Figure 48. Wait about 1 min before you inject again (step 13.b, below).
- 13. Perform the following repetitive sequence to obtain a total of four consecutive loop injections of reserpine.
- a. Select the Load option button, and overfill the injector loop with the 125 fg/ μ L solution of reserpine.
- b. Select the Inject option button to inject the reserpine solution into the ESI source, and then observe the Spectrum view.
- c. Wait 1 min before performing the next injection.
- d. Perform step 13.a through step 13.c three more times.
- 14. Click **Close** in the Divert/Inject Valve dialog box to return to the Tune Plus window.
- 15. Click **Stop** in the Acquire Data dialog box to end the data acquisition. Click **Cancel** to close the Acquire Data dialog box.

5 Acquiring ESI Sample Data Using the Tune Plus Window

Acquiring MS Data in the SIM Scan

19 C:\Xcalibur\methods\reserpine.LTQTune - Tune Plus File View Control ScanMode Display Setup Diagnostics H	eh	_ <u> 5 </u> ×
		LTO XL
	<u>ብ</u> ተ	
F: ITMS + p ESI SIM ms [609.20]		
100 96 90 85 80 76 70	Acquire Data Folder: C:VXcalbur\Data Folder: C:VXcalbur\Data File Name: reserptine C:CCalbur\Data File Name: reserptine File Name: reserve: r	? ×
60 60 4fsuatu etu etu etu etu etu etu etu etu etu e	Sample Name: reserpine Signs 10 Comment: SIM, ESI, 50pg loop Go Contact Closure Divgt Valve Start Mode OK Cancel Help	Peuse View Inst. Setup
30 26 20 16 10 5 0 606 607	609.7 608 609 610 m/z	613 614 615
For Help, press F1		NUM 6/11/2003 2:04 PM

Figure 48. Tune Plus window, showing the SIM spectrum of reserpine during analysis by loop injection

Review the mass spectrum and chromatogram in the Xcalibur Qual Browser window. See Figure 49.

For more information about reviewing the data you acquired using the LTQ XL MS detector with the Xcalibur data system, refer to the *Xcalibur Getting Productive: Qualitative Analysis* manual.

Acquiring MS Data in the SIM Scan



Figure 49. Qual Browser window, showing loop injections of reserpine in the Chromatogram view (left) and showing m/z 609 in the Spectrum view. Note that the injections occur at intervals of approximately 1 min

5 Acquiring ESI Sample Data Using the Tune Plus Window Acquiring MS Data in the SIM Scan

Chapter 6 Setting Up the Ion Source for Acquiring Data in APCI/MS/MS Mode

This chapter provides information on setting up the ion source for acquiring data in the APCI/MS/MS mode.

This chapter contains the following sections:

- Removing the ESI Probe
- Removing the Ion Max Ion Source Housing
- Removing the Ion Sweep Cone (optional)
- Installing the Corona Needle
- Installing the Ion Max Ion Source Housing
- Installing the APCI Probe

Removing the ESI Probe



Remove the ESI probe from the Ion Max ion source housing

- 1. Place the LC/MS system in Standby:
 - a. Stop the flow of solvent to the ESI source:
 - i. On the Control / Scan Mode toolbar, click **Inlet Direct Control** to display the Inlet Direct Control dialog box.
 - ii. In the Direct Control Panel group box, click the **Pump Off** button to stop the flow of solvent.
 - b. Click **On/Standby** on the Control / Scan Mode toolbar to place the MS detector in Standby.
- 2. Disconnect the sample transfer tubing from the stainless steel ZDV fitting (grounding union). See Figure 50.
- 3. Remove the 8 kV cable from the ESI needle high voltage receptacle shown in Figure 50 as follows:



CAUTION Handle the 8 kV cable with care.

- a. Unlock the cable by rotating the locking ring counter-clockwise.
- b. Unplug the 8 kV cable from the ESI needle high voltage receptacle.



Figure 50. Ion Max ion source housing with ESI probe installed

- 4. Disconnect the Auxiliary Gas fitting (green) from the auxiliary gas inlet (A) on the probe manifold. (Figure 50.)
- 5. Disconnect the Sheath Gas fitting (blue) from the sheath gas inlet (S) on the probe manifold.
- 6. Remove the stainless steel ZDV fitting (Grounding Union) from the grounding bar on the ion source housing.
- 7. Unlock the probe locking lever by rotating the lever open to its widest position.
- 8. Carefully pull the probe straight back in the port in the housing until it meets with the slot in the API interlock block. The guide pin on the probe manifold will prevent you from rotating the probe until the pin is aligned with the slot in the API interlock block. Once the probe is all the way back and aligned with the slot, turn the probe 45 degrees

counter-clockwise to free the probe from the alignment notch. Be careful not to break the fused-silica sample tube or PEEK safety sleeve.

- 9. Pull the probe straight out to remove it from the ion source housing.
- 10. Store the ESI probe in its original shipping container.

The ESI probe is properly removed from the Ion Max ion source housing.

Removing the Ion Max Ion Source Housing

You need to remove the ion source housing to access the ion sweep cone.

Note Disconnect any external liquid lines connected to the ion source housing before removing the ion source housing.

Remove the Ion Max ion source housing

- 1. Remove the drain tube from the ion source housing drain. See Figure 51.
- 2. Rotate the ion source housing locking levers 90 degrees to release the ion source housing from the ion source mount assembly.
- 3. Remove the ion source housing by pulling the housing straight off of the ion source mount assembly.
- 4. Place the housing in a safe location for temporary storage.

The Ion Max ion source housing is now properly removed. If you want to remove the ion sweep cone, go to the next section, Removing the Ion Sweep Cone. Otherwise, go to the topic "Installing the Corona Needle" on page 107.





Removing the lon Sweep Cone



Use of the ion sweep cone is optional. If you do not need to use the ion sweep cone it can be removed.

Remove the ion sweep cone

1. Wear a pair of talc-free gloves. Talc-free gloves are required so that the ion sweep cone is not contaminated.

CAUTION AVOID BURNS. At operating temperatures, the ion transfer capillary can severely burn you! The ion transfer tube typically operates between 200 and 400 °C. Always allow the ion sweep cone to cool to room temperature (for approximately 20 min) before you touch or remove this component. Always be careful not to touch the entrance end of the ion transfer capillary when it is exposed.

- 2. Grasp the outer ridges of the ion sweep cone and pull the cone straight off of the API cone seal.
- 3. Store the ion sweep cone in its original shipping container.

The ion sweep cone is now properly removed.

If you need to install the corona needle, go to the next section, Installing the Corona Needle. Otherwise, go to the section, "Installing the Ion Max Ion Source Housing" on page 108.

Installing the Corona Needle



Install the corona needle from the rear of the Ion Max ion source housing

CAUTION AVOID INJURY. The corona discharge needle is very sharp and can puncture your skin. Handle it with care.

1. Using needle nose pliers, grasp the corona needle and push the needle straight into the needle contact in the Ion Max ion source housing. Be careful not to scratch the needle. See Figure 52.





2. Make sure the tip of the needle is aligned with the path of travel between the APCI probe and the ion source interface on the instrument.

The corona needle is now properly installed.

Installing the Ion Max Ion Source Housing

Reinstall the Ion Max ion source housing

1. Carefully align the two guide pin holes on the rear of the ion source housing with the ion source housing guide pins on the MS detector, and carefully press the ion source housing onto the ion source mount. See Figure 53 and Figure 54.



Figure 53. Rear view of the Ion Max ion source housing



Figure 54. Ion source mount showing ion source housing guide pins

2. Rotate the ion source housing locking levers 90 degrees to lock the ion source housing onto the ion source mount assembly.

CAUTION Prevent solvent waste from backing up into the ion source housing and MS detector. Always ensure that liquid in the drain tube is able to drain to a waste container and that the outlet of the drain tube is above the level of liquid in the waste container.



CAUTION Do not vent the API source drain tube (or any vent tubing connected to the waste container) to the same fume exhaust system to which you have connected the forepumps. The analyzer optics can become contaminated if the API source drain tube and the (blue) forepump exhaust tubing are connected to the same fume exhaust system.

Your laboratory must be equipped with at least two fume exhaust systems. Route the (blue) forepump exhaust tubing to a dedicated fume exhaust system. Route the drain tube from the API source to a waste container. Vent the waste container to a dedicated fume exhaust system

3. Reinstall the ion source drain tube as follows:

- a. Connect the 1-in. ID Tygon tubing to the ion source housing drain fitting.
- b. Attach the free end of the hose to a waste container. Ideally, the waste container should be vented to a fume exhaust system.

The Ion Max ion source housing is now properly installed on the MS detector.

Installing the APCI Probe



Install the APCI probe into the Ion Max ion source housing

1. Connect the 8 kV cable to the corona needle high voltage receptacle as follows:

CAUTION Handle the 8 kV cable with care.

- a. Plug the 8 kV cable into the corona needle high voltage receptacle on the right side of the top of the ion source housing.
- b. Lock the cable by rotating the locking ring clockwise.
- 2. Be sure to unlock the probe locking lever (widest open position) before attempting to install the probe.
- 3. Insert the APCI probe into the port in the ion source housing, align the guide pin on the probe body at a 45 degree angle from the API interlock block. See Figure 55
- 4. Push the probe into the port until the guide pin meets with the probe collar on the ion source housing.
- 5. Turn the probe 45 degrees clockwise and align the guide pin with the slot in the API interlock block (you might need 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. Seat the probe all the way down into the alignment notch.



Figure 55. APCI probe

- 7. Lock the probe in place by rotating the probe locking lever towards the front of the housing; closing the probe locking lever towards the rear of the ion source housing might make it difficult to unlock.
- 8. Unplug the vaporizer heater cable from the ESI interlock plug on the ion source housing.
- 9. Connect the vaporizer heater cable to the vaporizer heater cable socket on the APCI probe.
- 10. Connect the sheath gas fitting (blue) to the sheath gas inlet (S) on the probe. (See Figure 55)
- 11. Connect the auxiliary gas fitting (green) to the auxiliary gas inlet (A) on the probe. (Figure 55)
- 12. Connect the sample transfer line to the APCI probe inlet.

The APCI probe is now properly installed in the Ion Max ion source housing.



CAUTION Prevent solvent waste from backing up into the ion source and MS detector. Always ensure that liquid in the drain tube is able to drain to a waste container and that the outlet of the drain tube is above the level of liquid in the waste container.

Leave the LC/MS system in Standby and go to Chapter 7: "Optimizing the MS Detector with Your Analyte in APCI/MS Mode".

Chapter 7 Optimizing the MS Detector with Your Analyte in APCI/MS Mode

This chapter provides information on optimizing the tune of your MS detector in the APCI/MS high flow mode. It is not necessary to recalibrate the MS detector when you switch to APCI/MS operation. You can use the calibration settings you obtained from the successful automatic calibration procedure you performed in the ESI/MS mode.

For APCI/MS operation you simply open a default Tune Method located in your C:\Xcalibur\methods folder, in this case APCIhighflow.LTQTune. From this starting point, you optimize automatically the tube lens voltage for your particular analyte. The capillary voltage and ion transfer capillary temperature may then be optimized manually to enhance ion transmission

Note The following procedures assume that you are familiar with your LTQ XL instrument and the Tune Plus window. For more information, refer to the LTQ XL online Help, *LTQ XL Getting Connected*, or the *LTQ XL Hardware Manual*.

Ensure that you have completed the procedures in the topics Tuning and Calibrating Automatically in the ESI/MS Mode and Setting Up to Acquire Data in the APCI/MS Mode.

This chapter includes the following sections:

- Setting Up the Inlet for Tuning Using High-Flow Infusion
- Setting Up the MS Detector for APCI/MS Operation
- Optimizing the Tune of the MS Detector Automatically in APCI/MS Mode
- Saving the APCI/MS Tune Method
- Cleaning the MS Detector after Tuning in APCI Mode

Setting Up the Inlet for Tuning Using High-Flow Infusion

Plumbing connections for APCI/MS sample introduction from the syringe pump into the LC solvent flow

1. Connect a 4 cm (1.5 in) segment of Teflon tubing with a (brown) fingertight fitting and a (brown) ferrule to the (black) LC union. See Figure 56.





- Load a clean, 500-μL Unimetrics syringe with 450 μL of a 125 fg/μL solution of reserpine or your analyte of interest. (See Appendix A: "Sample Formulations" for a procedure for making the reserpine tuning solution.)
- 3. Insert the needle of a syringe into the segment of Teflon tubing and place the syringe in the syringe holder of the syringe pump.
- 4. Connect a fused-silica infusion line from the (black) LC union to the (black) LC Tee union, as follows. See Figure 57.
 - a. Connect the infusion line with a (brown) fingertight fitting and a (brown) ferrule to the free end of the LC union.
 - b. Connect the other end of the infusion line with a (red) fingertight fitting and a (brown) ferrule to the side arm of the LC Tee union.



Figure 57. APCI/MS plumbing connections for the LC Tee union

Note To cut the PEEK tubing used to connect your LC to the divert/inject valve and the divert/inject valve to the APCI source, use a PEEK tubing cutter. This ensures that the tubing is cut straight. In addition, make sure your LC fittings, ferrules, and PEEK tubing are installed properly. By using these precautions, you prevent void (dead) volumes. The exclusion of void volumes is critical to microbore LC. Also, void volumes affect the quality of the MS detector signal.

- 5. Connect a segment of PEEK tubing from the (black) LC Tee union to the APCI LC inlet, as follows. See Figure 57.
 - a. Use a PEEK tubing cutter to cut a 4 cm (1.5 in.) length of the PEEK tubing.
 - b. Connect the PEEK tubing with a (brown) fingertight fitting and a (brown) ferrule to a free end of the (black) LC Tee union.
 - c. Connect the other end of the PEEK tubing with a (red) fingertight fitting and a (brown) ferrule to the LC inlet located on the APCI probe.

- 6. Connect an appropriate length of PEEK tubing (transfer line from the divert/inject valve) from the divert/inject valve to the LC Tee union, as follows. (See Figure 57.)
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and a (stainless steel) ferrule to port 3 of the divert/inject valve.
 - b. Connect the other end of the PEEK tubing with a (brown) fingertight fitting and a (brown) ferrule to the free end of the LC Tee union.
- 7. Connect an appropriate length of PEEK tubing (transfer line from the LC) from the divert/inject valve to the LC, as follows:
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and a (stainless steel) ferrule to port 2 of the divert/inject valve.
 - b. Connect the other end of the PEEK tubing with a proper fitting and a ferrule to the outlet of the LC.
- 8. Connect an appropriate length of PEEK tubing (waste line) from the divert/inject valve to a waste container, as follows:
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and a (stainless steel) ferrule to port 1 of the divert/inject valve.
 - b. Insert the other end of the PEEK tubing in a suitable waste container.

The LC plumbing connections are now properly made for APCI/MS sample introduction from the syringe pump into solvent flow from an LC.

Setting Up the MS Detector for APCI/MS Operation



Set up the MS detector for APCI/MS operation

- 1. In Tune Plus, click **On/Standby** to take the MS detector out of Standby mode and turn it On. The MS detector begins scanning, and the LTQ XL MS detector applies high voltage to the corona needle and shows a real-time display in the Spectrum view.
- 2. Open the *APCIhighflow.LTQTune* Tune Method, the Tune Method for high-flow APCI operation, as follows:
 - a. Choose **File > Open** to display the Open dialog box.
 - b. Scroll down until you see the folder *C:\Xcalibur\methods*. Then, select the file *APCIhighflow.LTQTune*.
 - c. Click **OK** to open the file. LTQ XL MS detector downloads the Tune Method parameters to the MS detector.
- 3. Verify that the LTQ XL MS detector opened the Tune Method, as follows:
 - a. On the Instrument Setup toolbar, click **API Source** to open the APCI Source dialog box. See Figure 58.

APCI Source		×
		Actual
⊻aporizer Temp (*C):	450.00	449.90
<u>S</u> heath Gas Flow Rate (arb):	80 ÷	78.91
Aux <u>G</u> as Flow Rate (arb):	20 ÷	18.96
S <u>w</u> eep Gas Flow Rate (arb):	0 +	0.00
Discharge Current (µA):	5.00	3.22
Discharge Voltage (kV) :		4.53
<u>C</u> apillary Temp (*C):	200.00	200.20
Capillary Voltage (V):	9.00	8.56
<u>T</u> ube Lens Offset (V):	100.00	100.15
Apply OK C	ancel <u>I</u>	<u>H</u> elp

Figure 58. APCI Source dialog box, showing the proper settings for a typical high flow experiment





- b. Verify that the settings in your dialog box are similar to those shown in Figure 58.
- c. Click **OK** to close the APCI Source dialog box.
- 4. Define the scan parameters for tuning the MS detector in the APCI/MS mode, as follows:
 - a. On the Control/Scan Mode toolbar, click **Define Scan** to open the Define Scan dialog box. Figure 59 (If your dialog box appears different from the one shown in the figure, it is probably because the advanced settings are not displayed. You can turn on the advanced settings as follows: In Tune Plus, choose **ScanMode**, and then click **Advanced Scan Features** to select the option.)
 - b. In the Scan Description group box, in the Mass Range list box, select **Normal** to allow for a selection of mass ranges between m/z 150 to 2000.
 - c. In the Scan Rate list box, select **Normal** to specify a normal scan rate.
 - d. In the Scan Type list box, select **SIM** to specify a selected ion monitoring scan.
 - e. In the Scan Time group box, in the Microscans spin box, type 1 to set the total number of microscans to 1.
 - f. In the Max. Inject Time spin box, type *200.000* to specify a 200 ms maximum injection time.

Define Scan										>
Scan History: ITMS + c SIM ms [608.20-610.20]										
Scan Description	EM	MSn Settings								
Mass Range: Normal 💌		n	Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)	#	Center Mass (m/z)	Width (m/z)
		2		1.0	20.0	0.250	10.000		609.20	2.00
Scan <u>Type:</u> SIM		3		1.0	20.0	0.250	10.000		2	
		4		1.0	20.0	0.250	10.000		3	
		5		1.0	20.0	0.250	10.000		1	
Scan Time		6		1.0	20.0	0.250	10.000		5	
Microscans: 1		7		1.0	20.0	0.250	10.000		6	
May Inject Time (me): 200,000		8		1.0	20.0	0.250	10.000		7	
		9		1.0	20.0	0.250	10.000		3	
	T	10		1.0	20.0	0.250	10.000		3	
Source Fragmentation					1	nput: Center	Width 💌			
Apply OK Cancel Help Injection RE Activation						Acti <u>v</u> ation				

Figure 59. Define Scan dialog box, showing typical settings for APCI/MS operation

- g. In the Source Fragmentation group box, confirm that the On check box is clear () to specify that the ion source fragmentation option is turned off.
- h. In the Scan Ranges group box, in the Input list box, select *Center/Width* to make available the Center Mass and Width text boxes in the Scan Ranges table.
- i. In the Scan Ranges group box, in the Scan Ranges table, in the Center Mass text box, type 609.20 to set the center mass for the scan range to m/z 609.20.
- j. In the Width text box, type 2.00 to set the width of the scan range to m/z 2.00.
- k. Verify that the settings in your Define Scan dialog box are the same as those shown in Figure 59.
- 1. Click **OK** to apply the MS detector scan parameters and to close the Define Scan dialog box.
- 5. On the Control/Scan Mode toolbar, click **Centroid/Profile** to toggle the data type to centroid. (The picture on the button should be the same as that shown here).

7 Optimizing the MS Detector with Your Analyte in APCI/MS Mode

Setting Up the MS Detector for APCI/MS Operation



6. Click the **Positive/Negative** button to toggle the ion polarity mode to positive. (The picture on the button should be the same as that shown here).

You have now completed setting up your MS detector for APCI/MS operation.

Optimizing the Tune of the MS Detector Automatically in APCI/MS Mode

You can optimize the tune of the MS detector automatically for APCI operation.

The most important parameters that affect the signal quality during APCI/MS operation are vaporizer temperature, ion transfer tube temperature, API gas flows, and solution flow rate. If any one of these parameters changes, you must re-optimize MS detector parameters. (You can use the Semi-Automatic tune procedure to tune the MS detector on individual parameters.)

Use the following procedure to automatically optimize the MS detector on the reserpine peak at m/z 609.2 at your particular flow rate, for example, 400 μ L/min. (See Table 3 on page 12 for guidelines about flow rates and temperatures.)

Automatically optimize the MS detector

- 1. On the Control/Scan Mode toolbar, click **Tune** to display the Automatic tuning page. See Figure 60.
- 2. In the What to Optimize On group box, select the Mass option button to make active the Mass spin box.
- 3. In the Mass spin box, type **609.2** to specify that you want to tune on the peak at m/z 609.2.
- 4. Verify that the Divert/Inject valve is in the Detector position, as follows:
 - a. Click the **Divert/Inject Valve** button to open the Divert/Inject Valve dialog box.
 - b. Select the **Detector** option button, and then click **Close** to return to Tune Plus.
- 5. Start the automatic tuning procedure from the Tune dialog box, as follows:
 - a. Click **Start**. A message box displays the following message:

Please ensure that the 500 microliter syringe is full.

Ensure the syringe pump contains at least 450 μL of the 125 fg/ μL reserpine tuning solution.

b. Click **OK** to close the message box, and return to the Tune Plus window.



7 Optimizing the MS Detector with Your Analyte in APCI/MS Mode

Optimizing the Tune of the MS Detector Automatically in APCI/MS Mode

П	
	f_/
	+

6. On the File/Display toolbar, click the Graph View button to display the view.

Tune
Automatic Semi-Automatic Manual Collision Energy
What to Optimize On
C Base Peak
• Mass (m/z): 609.20
Status
<u>S</u> tart Cancel <u>Print</u> <u>H</u> elp

Figure 60. Tune dialog box, showing the Automatic tuning page

7. Observe the Tune Plus window and the Tune dialog box. While automatic tuning is in progress, the LTQ XL MS detector displays various tests in the Spectrum and Graph views in the Tune Plus window and displays various messages in the Status group box in the Tune dialog box. Your Tune Plus window should now look similar to the one shown in Figure 61.

You have now successfully tuned the MS detector in APCI/MS mode for the compound reserpine (or your analyte of interest). Leave the LC pumps on (with a flow rate of approximately 400 μ L/min), and leave the Tune Plus window open with *APCIhighflow.LTQTune* to complete the next section, Saving the APCI/MS Tune Method.



Figure 61. Tune Plus window with the Tune dialog box, showing the Automatic tuning page

Saving the APCI/MS Tune Method

You can save the settings you just obtained in a Tune Method specific to your particular analyte and solvent flow rate. (In this case, you save settings obtained using reserpine.) You can recall the Tune Method and use it as a starting point for optimizing the MS detector on reserpine at a different flow rate.

Note Note. Save the Tune Method while the MS detector is still On.

Save the APCI/MS Tune Method

1. Choose File > Save As to display the Save As dialog box. See Figure 62.

Select the C:\Xcalibur\methods folder.

2. Click the **File Name** text box, and then enter **APCImyTune** to name the Tune Method APCImyTune.LTQTune.

Save As			?×
Save jn: 🗀	methods	- t	* III •
APCI.LTQ1 APCIHighF AutoTune.l APCILowFI APCITune. APCITune.l AutoTune.l	Tune Tow.LTQTune LTQTune low.LTQTune LTQTune LTQTune Tune	UESI.LTQTune ESIHighFlow.LTQTune ESILowFlow.LTQTune ESITune.LTQTune	
File <u>n</u> ame:	APCImyTune		<u>S</u> ave
Save as <u>type</u> :	Tune Files (*.LTQ	Tune)	Cancel
No file selec	mation		



3. Click **Save** to save the Tune Method and close the Save As dialog box. Note that the Tune Method is named *APCImyTune.LTQTune*.

Before you acquire data, go to the next section, Cleaning the MS Detector after Tuning in APCI Mode.

Cleaning the MS Detector after Tuning in APCI Mode





Use the following procedure to clean the MS detector after tuning on your analyte of interest.

1. Click **On/Standby** to put the MS detector in Standby mode. When the MS detector is in Standby, the LTQ XL MS detector turns off the vaporizer heater, corona discharge voltage, and syringe pump. The MS detector stops scanning, and the LTQ XL MS detector freezes the displays for the Spectrum and Graph views.

CAUTION Always place the MS detector in Standby (or Off) before you open the API source to atmospheric oxygen. The presence of oxygen in the ion source when the MS detector is On could be unsafe. (The LTQ XL MS detector automatically turns Off when you open the API source, however, it is best to take this added precaution.)

- 2. Remove the syringe from the syringe pump holder, as follows:
 - a. Squeeze the blue buttons, and pull back on the syringe pump handle to free the syringe.
 - b. Remove the syringe from the holder.
 - c. Disconnect the tip of the syringe needle from the Teflon tubing.
- 3. Clean the syringe thoroughly, as follows:
 - a. Clean the syringe with a solution of 5% formic acid in water.
 - b. Rinse the syringe with a solution of 50:50 methanol:water.
 - c. Use acetone to rinse the syringe. (Repeat this step several times.)



CAUTION AVOID BURNS. The APCI vaporizer heater can reach temperatures of 600 ×C. Always allow the APCI probe to cool to ambient temperature, for approximately 20 min, before handling or removing the APCI probe from the APCI flange.



CAUTION AVOID INJURY. The corona discharge needle is very sharp and can puncture your skin if you handle it without caution.

4. Remove the Ion Max ion source housing as described in the topic "Removing the Ion Max Ion Source Housing" on page 33. 5. Flush the sample transfer line, sample tube, and APCI probe thoroughly with a solution of 5% formic acid in water (or with another appropriate solvent), as follows:

Note The solvent that you use to flush the sample transfer line, sample tube, and APCI probe assembly depends on the solvent system you use to dissolve your samples. For example, if you are using a buffered solution of a high concentration, an acidic solution is appropriate.

- a. Fill a clean, 250 µL Unimetrics syringe with an appropriate solvent.
- b. While holding the plunger of the syringe in place, carefully insert the needle of the syringe into the free end of the Teflon tube.
- c. Flush the sample transfer line, sample tube, and APCI probe with the solution by slowly depressing the syringe plunger. Visually check that the solution is exiting the tip of the APCI probe on the inside of the probe assembly. Use a lint-free tissue to gently remove the excess solution as it exits the probe.
- d. Remove the needle of the syringe from the Teflon tube.
- 6. Repeat step 5 with a solution of 50:50 methanol:water.
- 7. Reinstall the Ion Max ion source housing as described in "Installing the Ion Max Ion Source Housing" on page 36.

If you plan to run analytical samples in APCI mode, go to Chapter 8: "Acquiring APCI Sample Data Using the Tune Plus Window".

Chapter 8 Acquiring APCI Sample Data Using the Tune Plus Window

This chapter provides information on acquiring LC/APCI/MS sample data using the Tune Plus window. This experiment uses reserpine but you can follow the same procedure with your analyte of interest.

Note The following procedures assume that you are familiar with your LTQ XL instrument and the Tune Plus window. If you need information, refer to the LTQ XL online Help, *LTQ XL Getting Connected*, and/or *LTQ XL Hardware Manual*. Verify that the procedures in Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode" and Chapter 7: "Optimizing the MS Detector with Your Analyte in APCI/MS Mode" have been completed .

This chapter contains the following topics:

- Setting Up to Introduce Sample by Loop Injection into Solvent Flow from an LC
- Acquiring APCI Data in the SIM Scan Mode

Setting Up to Introduce Sample by Loop Injection into Solvent Flow from an LC

This topic provides information on how to introduce sample by loop injection from an LC into solvent flow.

Plumbing connections

- 1. Connect an appropriate length of (red) PEEK tubing (transfer line from the divert/inject valve) from port 3 of the divert/inject valve to the (stainless steel) sample inlet fitting on the APCI probe. See Figure 63 and Figure 64.
- 2. Connect a 5 μ L sample loop with (stainless steel) set nuts and (stainless steel) ferrules to ports 1 and 4 of the divert/inject valve.
- 3. Connect an appropriate length of PEEK tubing (transfer line from the LC) from the divert/inject valve to the LC, as follows:
 - a. Connect a length of the PEEK tubing with a (stainless steel) nut and (stainless steel) ferrule to port 2 of the divert/inject valve.
 - b. Connect the other end of the PEEK tubing with a proper fitting and a ferrule to the outlet of the LC.



Figure 63. Divert/Inject valve, showing the correct setup for analysis by loop injection and showing the flow of liquid through the valve in the Load and Inject positions


Figure 64. APCI probe

- 4. Connect an appropriate length of PEEK tubing (waste line) from the divert/inject valve to a waste container, as follows:
 - a. Connect a length of PEEK tubing with a (stainless steel) nut and (stainless steel) ferrule to port 6 of the divert/inject valve.
 - b. Insert the other end of the PEEK tubing into a suitable waste container.

The MS detector is now set up to introduce sample by loop injection into solvent flow from an LC.

scan mode.

Acquiring APCI Data in the SIM Scan Mode





Note The LTQ XL MS detector automatically saves the acquired data on the hard disk.

Acquire a file of reserpine data in the selected ion monitoring (SIM)

- 1. If you have not already done so, in Tune Plus, click **On/Standby** to take the MS detector out of Standby mode and turn it On. The MS detector begins scanning, and the LTQ XL MS detector applies high voltage to the corona needle and shows a real-time display in the Spectrum view.
- 2. Ensure that the Centroid data type is selected. (The picture on the button should be the same as that shown here.)
- 3. Ensure that the scan parameters are defined to acquire SIM data for reserpine (or your analyte of interest), as follows:



- a. Click the **Define Scan** button to open the Define Scan dialog box. Figure 65
- b. Verify that the values in your dialog box are the same as those in Figure 8-3. Then, click **OK**.

Define Scan										
Scan <u>H</u> istory: ITMS + c SIM ms [60	08.20-61	10.20]								- 🖬 🛍
Scan Description	_ MSn	Settings						Sca	n Ranges —	
Mass Range: Normal 💌	n	Parent Mass (m/z)	lsolation Width (m/z)	Normalized Collision Energy	Activation Q	Activation Time (ms)		#	Center Mass (m/z)	Width (m/z)
	2		1.0	20.0	0.250	10.000		1	609.20	2.00
Scan <u>T</u> ype: SIM 🗾	3		1.0	20.0	0.250	10.000		2		
	4		1.0	20.0	0.250	10.000		3		
	5		1.0	20.0	0.250	10.000		4		
Scan Time	6		1.0	20.0	0.250	10.000		5		
Microscans: 1	7		1.0	20.0	0.250	10.000		6		
May Inject Time (ma): 200.000	8		1.0	20.0	0.250	10.000		7		
	9		1.0	20.0	0.250	10.000		8		
	10		1.0	20.0	0.250	10.000		9		
Source Fragmentation		⊻ideband Ac	tivation					10 <u>[</u> n	put: Center	Width 💌
	Apply	y 🗌 🗖	ОК	Cancel	Hel	p	Inj	jectio	n R <u>E</u> /	Activation

Figure 65. Define Scan dialog box, showing typical settings for acquiring data in the SIM scan mode

4. Turn on your LC pump, and specify an appropriate flow rate of $400 \ \mu$ L/min, for example. Verify that your system is free of leaks.

- 5. On the Control/Scan Mode toolbar, click **Acquire Data** to open the Acquire Data dialog box. See Figure 66.
- 6. Specify the acquisition parameters, as follows:
 - a. In the File Name box, type **reserpine** to specify a filename.
 - b. In the Sample Name box, type **reserpine** to specify the sample identity. If you are not using reserpine, type the name of your particular analyte.
 - c. In the Comment text box, type a comment about your experiment. For example, enter SIM, APCI, 50 pg, loop to specify the scan mode, ionization mode, sample amount, and/or method of sample introduction. The Xcalibur data system includes the comment on hard copies of your data.
 - d. In the Acquire Time box, select the Continuously option button to specify the continuous acquisition of data (until you stop the acquisition).
- 7. Leave the Acquire Data dialog box open during data acquisition, but move it to a corner of the Tune Plus window.
- 8. Click **Start** in the Acquire Data dialog box to begin acquiring data to the file reserpine3.raw. See Figure 67. The Acquisition Status group box displays the following message.

State: Acquiring Time (min):

Acquire Data			? ×
Fol <u>d</u> er: File Name:	C:\Xcalibur\Data	Acquire Time © Continuously	<u>S</u> tart
- Sample <u>N</u> ame:	reserpine	C Scans 10	Pause
<u>C</u> omment:	SIM, APCI, 50pg loop	<u>G</u> o to Standby when Finished	<u>V</u> iew
Instrument <u>M</u> ethod:	C:\Xcalibur\methods\MyMethod.meth	Acquisition Status State: Idle Time (min): 0.000	Inst. Setup
	Immediate Contact Closure Divert Valve OK Cancel Help		

Figure 66. Acquire Data dialog box, showing the acquisition status of the raw data file

8 Acquiring APCI Sample Data Using the Tune Plus Window

Acquiring APCI Data in the SIM Scan Mode

If C:\Xcalibur\methods\APCImyTune1TQTune-Tun File yew Control Scar@ode Display Setup Diagoes Image: Display Setup Diagoe	r Plus rs telo	LTO XL
Image: Second state of the second	Acquire Data Folger Folger Sample Name Comment Silm APCI, 50pg loop Bits Name Bits Name Comment Silm APCI, 50pg loop Bits Name Civicaliburing Statt Mode OK Cancel Help OK Cancel Help 608 609 600 610	21 ×1 Stop Pause Yiew Inst. Setup Inst. Setup 613
For Help, press F1	112	NUM 6/11/2003 2:04 PM

Figure 67. Tune Plus window, showing the SIM spectrum of reserpine during analysis by loop injection



- 9. Inject the reserpine solution into the APCI source from the Instrument Setup toolbar, as follows:
 - a. Click **Divert/Inject Valve** to display the Divert/Inject Valve dialog box. See Figure 68.

Divert/Inject V	alve	×
Load	Inject	
۲	0	
Detector	Waste	
Close	<u>H</u> elp	
		-

Figure 68. Divert/Inject Valve dialog box

- b. Select the **Load** option button, and overfill the 5- μ L injector loop with the 125 fg/ μ L solution of reserpine (or a solution of your analyte of interest).
- c. Select the Inject option button.
- 10. Observe the reserpine peak (m/z 609.2), or that of your analyte of interest, in the Spectrum view.
- 11. Perform the following repetitive sequence to obtain a total of four consecutive loop injections of reserpine in the SIM scan mode. Wait about 1 min between injections.
 - a. Select the Load option button to put the divert/inject valve in the Load position. Overfill the injector loop with the 125 fg/ μ L solution of reserpine.
 - b. Select the Inject option button to inject the reserpine solution into the APCI source. Then, observe the Spectrum view.
 - c. Wait 1 min before the next injection.
 - d. Repeat step 11.a through step 11.c three times
- 12. Click **Stop** in the Acquire Data dialog box to end the data acquisition.

Review the mass spectrum and chromatogram in the raw file you just acquired using the Xcalibur Qual Browser window. See Figure 69.

For more information about reviewing the data you acquire using the LTQ XL MS detector with the Xcalibur data system, refer to the *Xcalibur Getting Productive: Qualitative Analysis* manual.

Note If you want to acquire MS/MS Full scan data in APCI mode, see "Setting Up to Acquire MS/MS Data in the Full Scan Type" on page 84 for information about setting up the LTQ XL MS detector.



Figure 69. Qual Browser window, showing loop injections of reserpine in the Chromatogram view (left) and *m/z* 609 in the Spectrum view

Appendix A Sample Formulations

This appendix provides instructions for the preparation of several stock solutions. These solutions are used for tuning, calibrating, and demonstrating applications of the APCI / ESI system. Formulations for sample solutions in this appendix include:

- Caffeine, MRFA, and Ultramark 1621 Stock Solutions
- ESI Calibration Solution: Caffeine, MRFA, Ultramark 1621
- Reserpine

The caffeine, MRFA,Ultramark 1621, and Resperine needed to make the solutions are supplied with your chemical accessory kit. You can order replacement chemical accessory kits from Thermo Electron (Thermo Electron part #97000-62042).







CAUTION Store and handle all chemicals in accordance with standard safety procedures. The Material Safety Data Sheet (MSDS) describing the chemicals being used are to be freely available to lab personnel for them to examine at any time. Material Safety Data Sheets (MSDSs) provide summarized information on the hazard and toxicity of specific chemical compounds. MSDSs also provide information on the proper handling of compounds, first aid for accidental exposure, and procedures for cleaning spills or dealing with leaks. Producers and suppliers of chemical compounds are required by law to provide their customers with the most current health and safety information in the form of an MSDS. Read the MSDS for each chemical you use. Dispose of all laboratory reagents in the appropriate way (see the MSDS).

Potentially hazardous chemicals used in procedures throughout this manual include:

- Acetic acid
- Acetonitrile
- Methanol
- Reserpine
- Formic Acid

Caffeine, MRFA, and Ultramark 1621 Stock Solutions

For tuning and calibrating the ESI system, you use a solution of caffeine, MRFA, and Ultramark 1621 in an acetonitrile:methanol:water solution containing 1% acetic acid. You prepare the calibration solution from each of the following:

- Caffeine stock solution
- MRFA stock solution
- Ultramark 1621 stock solution

Note Vials of caffeine, MRFA, and Ultramark 1621 are included in the API accessory kit. These compounds are available from Sigma Chemical Co. The Sigma caffeine Product Number is C6035 (Caffeine Solution, drug standard 1mg/mL±5% in methanol). The Sigma MRFA Product Number is M1170 (MRFA=Met-Arg-Phe-Ala acetate salt). To order more of these compounds from Sigma, write or call:

Sigma Chemical Company P. O. Box 14508 St. Louis, Missouri, USA 63178-9916 (800) 325-3010 (in the USA) (905) 829-9500 (in Canada) (314) 771-3750 (outside the USA or Canada) www.sigmaaldrich.com **Note** A vial of Ultramark 1621 (neat liquid) is included in the API accessory kit. This compound is available from Lancaster Synthesis. The The structure of Ultramark 1621 is ahown below (x is 1, 2, or3).



The Lancaster Ultramark 1621 Catalog Number is 16698 (Ultramark 1621, Mass Spec Std.). To order more of this compound from Lancaster Synthesis, write or call:

Lancaster Synthesis, Inc. P.O. Box 1000 Windham, NH USA 03087-9977 (603) 889-3306, (800)-238-2324 (in the USA & Canada) +44 (0)1524 36101 (UK & International) www.lancastersynthesis.com



CAUTION AVOID EXPOSURE TO POTENTIALLY HARMFUL MATERIALS. Always wear protective gloves when you use solvents or corrosives. Refer to your supplier's Material Safety Data Sheet (MSDS) for the proper handling of a particular solvent. Contain waste streams, and use proper ventilation.



CAUTION AVOID EXPOSURE TO POTENTIALLY HARMFUL MATERIALS. Always wear safety glasses when you use solvents or corrosives. Refer to your supplier's Material Safety Data Sheet (MSDS) for the proper handling of a particular solvent. Contain waste streams, and use proper ventilation.

Stock Solution: Caffeine

A 1 mg/mL stock solution of caffeine in 100% methanol is provided with your LTQ XL system. This is also the concentration of the Sigma replacement mentioned in the Note above (Sigma #C6035).

Stock Solution: MRFA	Prepare a 1.5 mL stock solution of 166.7 pmol/ μL MRFA in 50:50 methanol:water
	 Obtain the vial of L-methionyl-arginyl-phenylalanyl-alanine acetate•H2O (MRFA) in your accessory kit. In this form, the MRFA sample has an average molecular weight of 523.7 u. Carefully weigh 2.6 mg of the MRFA sample.
	2. Dissolve the MRFA sample in a total volume of 1.0 mL of 50:50 methanol:water.
	3. Mix the solution (5.0 nmol/ μ L) thoroughly.
	4. Transfer 50 μL of the 5 nmol/μL solution into a clean polypropylene tube.
	5. Add 1.45 mL of 50:50 methanol:water to the tube.
	6. Mix this solution (166.7 pmol/µL) thoroughly.
	7. Label the tube MRFA stock solution.
Stock Solution: Ultramark	Prepare a 10 mL stock solution of 0.1% Ultramark 1621 in acetonitrile
1621	1. Obtain the vial of Ultramark 1621 in your accessory kit.
	2. Using a gas-tight syringe, measure out 10 μ L of Ultramark 1621. Special care in taking the sample may be needed as the Ultramark 1621 has a high viscosity. Dissolve the 10 μ L of Ultramark 1621 in 10 mL of acetonitrile.

- 3. Mix the solution thoroughly.
- 4. Label the vial Ultramark 1621 stock solution.

ESI Calibration Solution: Caffeine, MRFA, Ultramark 1621

Prepare 10 mL of ESI calibration solution

- 1. Pipet 200 μL of the stock solution of caffeine into a clean, dry 10 mL volumetric flask.
- 2. Pipet 100 μL of the stock solution of MRFA into the flask.
- 3. Pipet 100 μL of the stock solution of Ultramark 1621 into the flask.

Note Use only glass pipets or stainless steel syringes when measuring glacial acetic acid. Using plastic pipet tips causes contamination of acid stock solutions that can introduce contaminants in the calibration solution.

- 4. Pipet 100 μ L of glacial acetic acid into the flask.
- 5. Pipet 5mL of acetonitrile into the flask.
- 6. Bring the volume of the solution up to the 10 mL-mark on the flask with 50:50 methanol:water.
- 7. Mix the calibration solution thoroughly.
- 8. Transfer the solution to a clean, dry vial.
- 9. Label the vial *ESI Calibration Solution* and store it in a refrigerator until it is needed.

Reservine Use the following directions to prepare a stock solution of reserpine. Use serial dilutions of this stock solution to make a sample solution. **Stock Solution: Reserpine** Prepare a stock solution of 1 μ g/ μ L reserptine in 1% acetic acid-methanol 1. Obtain the 1 gram vial of reserpine in your accessory kit. (The average molecular weight of reserpine is 608.7 u). Weigh out 10 mg of reserpine and transfer the sample to a 10 mL volumetric flask. 2. Dilute the reserpine up to the 10 mL mark with a solvent of 1% acetic acid in methanol. 3. Ensure that the sample is thoroughly dissolved in solution. 4. Transfer the solution to a clean and dry light resistant vial. Label this vial as *Reserpine Stock Solution (1 mg/mL=1.64nmol/mL)*. It is best to use the Reserpine Stock solution as soon as it is made. If it is necessary to store the solution, keep it in a refrigerator until it is needed. **Reserpine Tuning Solution** Prepare 1 mL of either the Reservine Tuning Solution of 10 pg/ μ L (16.4 fmol/µL, step 9) or the Reserpine APCI Sample Solution of and Reserpine APCI 250 fg/ μ L (411 amol/ μ L, step 16) in 1% acetic acid-50:50 methanol:water. Sample Solution Prepare the appropriate Reserpine Solution 1. Pipet 100 μ L of the Stock Solution (1 μ g/ μ L) of reserpine into a clean polypropylene microcentrifuge tube. 2. Add 900 μ L of 1% acetic acid in 50:50 methanol:water to the tube. 3. Mix this solution (100 ng/ μ L=164 pmol/ μ L) thoroughly. 4. Transfer 10 μ L of the 100 ng/ μ L solution into a clean polypropylene tube. 5. Add 990 μ L of 1% acetic acid in 50:50 methanol:water to the tube. 6. Mix this solution (1 ng/ μ L=1.64pmol/ μ L) thoroughly. 7. Transfer 10 μ L of the 1 ng/ μ L solution into a clean polypropylene tube. 8. Add 990 μ L of 1% acetic acid in 50:50 methanol:water to the tube.

- 9. Mix this solution (10 pg/ μ L=16.4 fmol/ μ L) thoroughly. This is the final dilution for the Reserpine Tuning Solution. If you are stopping at this dilution, label it as *Reserpine Tuning Solution (10 pg/\muL)*. It best to use this solution immediately after it is made. Keep it in a light resistant container in a refrigerator if it must be stored. If you are making a Reserpine APCI Sample Solution, continue with step 10.
- 10. Transfer 100 μL of the 10 pg/ μL solution into a clean polypropylene tube.
- 11. Add 900 μL of 1% acetic acid in 50:50 methanol:water to the tube.
- 12. Mix this solution (1 pg/ μ L=1.64 fmol/ μ L) thoroughly.
- 13. Transfer 100 μL of the 1 pg/ μL solution into a clean polypropylene tube.
- 14. Add 300 μ L of 1% acetic acid in 50:50 methanol:water to the tube.
- 15. Mix this solution (250 fg/ μ L=411 amol/ μ L) thoroughly.
- 16. Label the tube *Reserpine APCI Sample Solution (250 fg/\muL)*. It is best to use this solution immediately after it is made. Keep it in a light-resistant container in a refrigerator if it must be stored.

Appendix B LTO XL High Mass Range Calibration

A high mass range calibration is necessary if you want to use your LTQ XL to analyze masses in the 2000 u-4000 u range. This appendix provides instructions for the high mass range calibration of the LTQ XL. The presentation in this Appendix assumes that you are familiar with the contents of Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode" and Appendix A: "Sample Formulations".

This Appendix Contains the following topics:

- High Mass Range Calibration Solution
- Normal Mass Range Calibration
- Enter Normal Mass Range Data into Tune Plus
- Tune on m/z 524.3
- High Mass Range Calibration Procedure
- Notes

High Mass Range Calibration Solution

The high mass range calibration solution is 70 ng/ μ L polypropylene glycol (PPG) in a solvent of 65/35 methanol/10mM sodium acetate. The PPG used for the calibration procedure is Aldrich #202347 as specified in the following Note.

Tip One possible formulation for this solution is 3.5mg of PPG diluted to 50mL with the 65/35 methanol/10mM sodium acetate solvent.

Note The procedures that follow assume that the high mass range calibrant is polypropylene glycol, number average molecular weight about 2700 (Mn~2700), Aldrich product number 202347. To order this compound from Sigma-Aldrich, write or call:

Sigma Chemical Company P. O. Box 14508 St. Louis, Missouri, USA 63178-9916 (800) 325-3010 (in the USA) (905) 829-9500 (in Canada) (314) 771-3750 (outside the USA or Canada) www.sigmaaldrich.com







CAUTION Store and handle all chemicals in accordance with standard safety procedures. The Material Safety Data Sheets (MSDS) describing the chemicals being used are to be freely available to lab personnel for them to examine at any time. Material Safety Data Sheets (MSDS) provide summarized information on the hazard and toxicity of specific chemical compounds. MSDSs also provide information on the proper handling of compounds, first aid for accidental exposure, and procedures for the remedy of spills or leaks. Producers and suppliers of chemical compounds are required by law to provide their customers with the most current health and safety information in the form of an MSDS. Read the material safety data sheets for each chemical you use.

Normal Mass Range Calibration

Before calibrating for the high mass range you must calibrate for the normal mass range. Use the calibration solution (ESI calibration solution) described in Appendix A: "Sample Formulations" and the procedures described in Chapter 3: "Tuning and Calibrating Automatically in the ESI/MS Mode" to perform the normal mass range calibration. The calibration solution contains caffeine as the low molecular mass component, MRFA as the mid-molecular mass component, and Ultramark 1621 as the high molecular mass component. Four pieces of information from this calibration are needed for the high mass range calibration:

- 1. The expected low mass (caffeine component=195.10)
- 2. The expected high mass (1822.00 peak of the Ultramark 1621 component)
- 3. Observed low mass (observed mass of the caffeine component)
- 4. Observed high mass (observed mass of the 1822 peak of the Ultramark 1621 component).

Enter Normal Mass Range Data into Tune Plus

Enter Normal Mass Range Data into Tune Plus

- 1. After carrying out the normal mass range calibration open the Tune Plus window (**Instrument Set Up>LXQ>Start Tune Plus**).
- 2. In the Tune Plus window, open the Diagnostics dialog box (**Diagnostics>Diagnostics**) shown in Figure 70.

Diagnostics	
ToolsTestsPlot readbackSet deviceRF tuneDevice calibrationDisplay settingsTogglesTriggersMass calibrationSystem evaluation	Manual coarse calibration Calibrate current scan type Mass list: User Expected low mass: 195.10 Expected high mass: 195.10 Dbserved high mass: 1819.30 Set AGC mode Execute Estimate other calibration modes using current full scan Execute Save Save
	OK Cancel Print Help

Figure 70. Tune Plus Diagnostics dialog box, normal mass range data

- 3. Click **Tools** in the upper left of the window. Click on Mass Calibration in the pane below the **Tools** button as shown in Figure 70.
- 4. Items mentioned in step 5 through step 7 below are inside of the Manual Coarse Calibration group box.
- 5. In the Calibrate Current Scan Type group box, choose User in the Mass list box. In the Expected Low Mass box type 195.10 and type 1822.00

in the Expected High Mass box. Type the Observed Low Mass and the Observed High Mass values in their respective text boxes. These latter values are obtained as discussed in the preceeding section: Normal Mass Range Calibration. After entering the observed mass values click the **Execute** button.

- 6. In the Estimate Other Calibration Modes Using Current Full Scan box, click **Execute**.
- 7. Click **Save** at the bottom right of the Manual Coarse Calibration window.
- 8. Click **OK** at the bottom of the Diagnostics dialog box.

Tune on m/z 524.3



Optimize the MS detector Tune on *m/z* 524.3u.

- 1. On the Control/Scan Mode toolbar, click on the Tune button to display the Tune dialog box.
- 2. If necessary, click on the Automatic tab to display the Automatic tuning page. See Figure 71.
- 3. In the What to Optimize On group box, select the Mass option button to make active the Mass spin box.
- 4. In the Mass spin box, type 524.3 to specify that the LXQ MS detector uses the peak at m/z 524.3 to optimize your tune.

Tune Tune				
Automatic Se	mi-Automatic M	anual Collision En	ergy	
What to Opt	imize On			
⊂ <u>B</u> ase P	eak n/z): 524.3	•		
Status				
	<u>S</u> tart	Cancel	Print	<u>H</u> elp



5. Click **Start**. A message box displays the following message:

Ensure that the 500 microliter syringe is full.

Ensure that the syringe pump contains at least 450 μ L of the ESI calibration solution as mentioned in the topic "Normal Mass Range Calibration" on page 149.

6. Click **OK** to close the message box, and return to the Tune Plus window.



- 7. In the File/Display toolbar, click the Graph View button to display the Graph view.
- 8. Observe the Tune Plus window and the Tune dialog box. While automatic tuning is in progress, the LXQ MS detector displays various tests in the Spectrum and Graph views in the Tune Plus window and displays various messages in the Status group box in the Tune dialog box.

High Mass Range Calibration Procedure

Calibrate the high mass range automatically in ESI mode

- 1. Load the syringe with PPG2700 solution.
- 2. From the Tune page go to Define Scan dialog box (see Figure 32 on page 76). In the Scan Description group box: a) choose **High** in the Mass Range spin box, b) choose **Normal** in the Scan Rate spin box, and c) choose **Full** in the Scan Type spin box. In the Scan Ranges box, type 2000 for the First Mass and 4000 for the Last Mass. Click **Apply**. The spectrum should look like that shown in Figure 72.
- 3. Do a two-point mass calibration. Use the Tune Plus Diagnostics dialog box (Figure 73). Access the Diagnostics dialog box as instructed in "Enter Normal Mass Range Data into Tune Plus" on page 150. Type the high mass values as shown in Figure 73. Click **Execute**, click **Save**, and close the window.
- 4. In the Tune Plus window, choose **Control>Calibrate** to open the Calibrate dialog box (Figure 74).



Figure 72. PPG2700 mass spectrum

Diagnostics	
ToolsTestsPlot readbackSet deviceRF tuneDevice calibrationDisplay settingsTogglesTriggersMass calibrationSystem evaluation	Manual coarse calibration Calibrate current scan type Mass list: User Expected Jow mass: 2654.21 Expected Jow mass: 2654.21 Expected high mass: 3700.06 Observed high mass: 3638 Set AGC mode Execute Estimate other calibration modes using current full scan Execute Save Save
	OK Cancel <u>Print</u> Help

Figure 73. Tune Plus Diagnostics dialog box, two point high mass calibration

Calibrate		×
High Mass Range Calibration	Mass Range: 🔿 Normal 🗭 High	
What to do	Calibration Mass List <u>Name:</u> PPG 2700 (factory)	
 Check Calibration What to cal/check ✓ Mass for Normal ✓ Mass for Iurbo ✓ Mass for Zoom ✓ Isolation Waveforms 	# Monoisotopic m/z Average m/z 1 2652.880000 2654.610000 2 3001.140000 3003.100000 3 3349.390000 3351.580000 4 3697.640000 3700.060000	
Status	Sa <u>v</u> e Save <u>A</u> s <u>D</u> elete	
Set instrument to Standby Wh	Cancel <u>Print</u> <u>H</u> elp	

Figure 74. Calibrate dialog box

- 5. For Mass Range, select the High button (upper right of the Calibrate dialog box). The High Mass Range Calibration tabbed page appears (as shown in Figure 74).
- 6. In the What To Do box, select the Calibrate button. Type a calibration mass list name in the Name list box inside the Calibration Mass List group box.

- If you want to use a previously saved mass list Name in the Calibration Mass List group box go to step 7.
- If you want to use a new mass list Name in the Calibration Mass List group box go to step 8.
- 7. To use a previously saved calibration mass list Name in the Calibration Mass List group box,
 - a. Select the Name list from the name of a previously saved calibration mass list. The calibration masses saved in the selected list appear in the table below the Name list box (in the Monoisotropic m/z and the Average m/z columns). The example in Figure 74 shows the calibration masses saved for the calibration mass list: PPG 2700 (factory).
 - b. Go to step 9.
- 8. To assign a new calibration mass list name,
 - a. You can use the factory default PPG2700 name that is in the Name list box or you may assign a new name for this mass list.
 - b. To assign a new name click on the **Save As** button. The Unique Mass List Name dialog box appears.
 - c. Type the new name in the Unique Mass List Name dialog box and click **OK** to accept the name. The dialog box closes and the new name appears in the calibration mass list Name list box.
 - d. Do not change the masses in the Calibration Mass List table.
- 9. In the 'What to cal/check' area, click the Mass for Normal, Mass for Turbo, Mass for Zoom, and the Isolation Waveforms check boxes as shown in Figure 74.
- 10. Specify the instrument state when the calibration is completed. Use the Set Instrument to Standby When Finished check box at the bottom left of the Calibrate dialog box.
 - If you want the LTQ XL to switch to Standby mode after the calibration, click the check box so that a check mark is present in the box.
 - If you want the LTQ XL to remain in the On mode after the calibration, clear the check box.
- 11. Start the calibration procedure.

- a. At the bottom of the Calibrate dialog box, click **Start**. A message box appears with the message: Please ensure that the 500 microliter syringe is full.
- b. Verify that the syringe contains at least 420μ L of the high mass range calibration solution.
- c. Click OK to close the message box and return to the Calibrate dialog box. Start the calibration.
- 12. Observe the Tune Plus window and the Calibrate dialog box. While the automatic calibration is in progress the Spectrum and Graph views display a variety of test results and the Status pane in the Calibrate dialog box displays status messages. The system will automatically use the average mass values for the Turbo Scan calibration. Monoisotopic masses will automatically be used for the normal, Zoom Scan and isolation waveform calibrations.
- 13. Observe the PPG2700 calibrations are complete. it may be necessary to repeat step 3-the two point mass calibration.

When all calibration items are completed, your LTQ XLis properly tuned and calibrated in the high mass range. At the conclusion of the procedure, you must clean the syringe and API source (change the tubing and syringe, clean the API source with acetone to remove the polypropylene glycol).

Notes

- 1. Two-point mass calibration is available for all high mass range modes accessed through the Estimate Other Calibration Modes Using Current Full Scan group box (in the Diagnostics dialog box, Figure 70).
- 2. Use the **Execute** button in the Estimate Other Calibration Modes Using Current Full Scan box to execute estimation for all other calibrations from the normal mass range-normal scan rate mass calibration. The other modes for which the calibrations get estimated include Zoom Scan, Turbo Scan, Enhanced Scan, Isolation Waveform, and High Mass Isolation Waveform calibrations.

This button changes when the High Mass Scan mode is selected. In this case, clicking on the **Execute** button causes the High Mass Isolation to be estimated from the Current Normal Mass Range Isolation calibration (assuming that it has been calibrated). This feature has been added to extrapolate the High Mass Isolation calibration for Initial calibration. As noted in item #1, a two point mass calibration is available for all other scan modes . This makes it possible to get sufficient convergence of various mass values to allow automatic calibration.

3. In order to calibrate with High Mass Range Zoom Scan, it may be necessary to reduce the Zoom Scan target from 3000 u to 1000u.

Index

A

Accept Optimized Value dialog box 91 Acquire Data dialog box displaying 95, 135 starting acquisition from 95, 135 stopping acquisition from 97, 137 acquiring data prerequisites for, in MS/MS mode 84 using Tune Plus 83, 131 APCI probe installing 111 APCI Source dialog box displaying 119 APCI vaporizer avoiding burns from (CAUTION) 29, 30, 66 disabling 40 APCI/MS setting up Finnigan LTQ for 119 using, discussion 5 APCI/MS operation, setting up 119 AutoTune file 49

B

Breadth Focus, in Ion Tree experiment, discussion 24 buffers, discussion 10

C

caffeine preparing calibration solution from 143 reordering (note) 140 spectral ion from 53 vial of stock solution description 141, 141 Caffeine Stock solution 141 Calibrate dialog box description 13 displaying 62 figure 64 messages in Status box 63 calibration discussion 13 ions 53 LTQ XL setup 47

procedure 62 solutions caution 74 preparing 143 syringe pump setup for 45 calibration solutions, preparation 143 CAUTIONS avoiding burns from APCI vaporizer 29, 30, 66, 128 from ion sweep cone 106, 106 from ion transfer capillary 35, 66 avoiding injury from corona discharge needle 31, 107 Cautions avoiding contamination of system 14 calibration solution, flow rate limit of 74 solvent waste backing up into ion source housing 37, 109, 113 venting API source 38, 109 cleaning Finnigan LTQ 65, 128 collision energy optimizing 89 setting, in Define Scan dialog box (note) 84 contamination caution 14 note 45 of solvents 143 corona discharge needle avoiding injury 31, 107 figure 32 installing 107 removing 31

D

Data-Dependent Ion Tree description 25 figure 24 Define Scan dialog box displaying 50, 75, 94, 120, 134 Depth Focus, discussion 24 dialog boxes Accept Optimized Value 91 APCI Source 119 Tune 124

E

ESI Calibration Solution Caffeine, MRFA, Ultramark 1621 143 ESI probe installing 39 ESI Source dialog box displaying 49, 53 figure 49, 59 viewing parameters 57 ESI/MS discussion 4 operational guidelines (table) 11 optimizing Tune Method 69 preparing calibration solution for 143 setting up Finnigan LTQ for 74 testing Finnigan LTQ 52 ESI/MS operation, setting up 74 experiment types data-dependent 17 discussion 16 General MS 16 Ion Mapping 21

F

flow rates discussion 11

G

Graph view displaying 78, 124, 153 freezing 65, 128 observing 63 guidelines LC/APCI/MS operational (table) 12 LC/ESI/MS operational (table) 12

Instrument Setup window description 16 displaying 28, 47 introducing sample 8 introducing samples into Finnigan LTQ, discussion 8 Ion Max ion source discussion 4 installing 108 removing 33, 105 Ion Optics dialog box

displaying 57 ion sweep cone avoiding burns from (CAUTION) 106, 106 description 34 installing 34 removing 106 ion transfer capillary avoiding burns from (CAUTION) 35, 66, 106 cleaning 66 discussion 11, 70 flushing 67 gaining access to 65 optimizing temperature of discussion 115, 123 note 53, 79 parameters affecting signal quality (note) 14 reinstalling 66 removing 66 sealing 66 temperature guidelines APCI operation 12 ESI operation 12 unsealing 67 Ion Tree experiment Breadth Focus 24 Depth Focus 24 isolation width note 87 optimizing 84 optimizing (note) 87

L

loop injection setup for 92, 132 LTQ LX cleaning for normal operation 128 LTQ XL calibrating 62 cleaning for normal operation 65 discussion 2 ESI/MS mode, testing 52 flow rates 11 introducing sample 8 operational guidelines 11 optimizing Tune Method, LC/ESI/MS operation 69 placing in Standby 28 setting up for APCI/MS operation 119 setting up for ESI/MS operation 74

Μ

MRFA preparing calibration solution from 143 reordering (note) 140 spectral ion from 53 MRFA Stock solution 142 MS/MS acquiring, data 84 discussion 3

Ν

nitrogen atomic, affect on ion current 6 sweep gas discussion 7 ensure flow to API source (CAUTION) 47 starting flow of 48 Normalized Collision Energy parameter 84

0

operational guidelines 11 optimizing auxiliary gas 7 collision energy 89 automatically 88, 89 discussion 84 note 84 saving value 90 Finnigan LTQ automatically 47 for LC/ESI/MS operation 69 ion transfer capillary temperature 56 isolation width 84 discussion 84 note 87 performance of data acquisition, discussion 13 pH of solvent 5 sweep gas 7 Tune Method automatically 77 discussion 57, 69 Tune Method (note) 57, 57 optimizing tune method, LC/ESI/MS operation 69

Ρ

PEEK tubing, cutting (note) 117 placing in Standby 28 procedures acquiring data using Tune Plus 83, 131 calibrating LTQ XL automatically 62 cleaning Finnigan LTQ after tuning and calibrating 128 cleaning LTQ XL after tuning and calibrating 65 installing APCI probe 111 corona discharge needle 107 ESI probe 39 Ion Max ion source 108 ion sweep cone 34 optimizing APCI tune 123 collision energy 89 ESI tune 77 isolation width 84 preparing calibration solutions 143 Ultramark 1621 stock solution 142 removing APCI probe 29 ESI probe 102 ion sweep cone 106 saving Tune Method 60, 81, 126 setting up hardware for tuning and calibrating 27 loop injection 92, 132 LTQ XL analyzing your analyte 74 LTQ XL for APCI/MS operation 119 syringe pump 45, 71 stopping LC flow 102 testing LTQ XL in ESI/MS mode 52 tuning and calibrating Finnigan LTQ 43 setting up for 47 tuning LTQ XL automatically in APCI/MS mode 123 in ESI/MS mode 56 using your analyte 69

0

quantitation experiments 16

R

Reserpine 144 reserpine acquiring, data, loop injection 83, 131 preparing solutions of 144 Reserpine APCI Sample Solution 144 Reserpine Stock solution 144 Reserpine Tuning Solution 144

S

Save As dialog box displaying 60, 81, 126 solutions, preparing calibration 143 Standby, placing Finnigan LTQ in 28 syringe pump description 8 installing syringe into 45, 71, 116 locating 45 plumbing 71, 116, 118 removing syringe from 65, 128 setting up for tuning and calibrating 45 starting 52, 57, 78, 85, 123, 153 stopping 65, 90, 128 using 13, 14 Syringe Pump dialog box displaying 86

Т

Tables ESI operational guidelines 11 **Tune Methods** opening predefined 49, 74, 119 Tune Plus window opening 28, 47, 74 placing Finnigan LTQ in Standby from 28 tuning results (figure) 124 using, to acquire data 83, 131 tuning automatically 56 discussion 13 ESI tune files (note) 49 LTQ XL for LC/ESI/MS operation 69 setup for 47, 74 LTQ XL (note) 57 syringe pump, setup for 45 voltages 56

U

Ultramark 1621 cleaning spray shield of 67 contamination from (in caution note) 14, 74 preparing calibration solution from 143 preparing stock solution of 142 reordering 140, 141 spectral ions from 53 Ultramark 1621Stock solution 142

V

voltages automatic optimization, description 43, 69

W

Wideband Activation discussion 3

X

Xcalibur displaying Home Page 28, 47, 47, 74 displaying Roadmap view 74

Ζ

ZoomScan, discussion 4, 20