

Notices

Manual Part Number

D0015483
June 2022

Copyright

© Agilent Technologies, Inc. 2022

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051

Software Revision

This guide is valid for the LTS 01.11 or later revision of the Agilent OpenLab ChemStation software for the Agilent InfinityLab LC/MSD Series and 6100 Series LC/MS, until superseded.

Software Manufacturing



Manufactured for
Agilent Technologies
5301 Stevens Creek Blvd
Santa Clara, CA 95051

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

In this Guide...

This guide presents a series of exercises to help you learn the basic operation of the Agilent OpenLab ChemStation software with your InfinityLab LC/MSD Series or 6100 Series LC/MS system.

1 Prepare for the Analysis

Use these exercises to prepare the LC, to dilute a sulfa demonstration sample, and to check the tune on the MS.

2 Set Up and Run a Scan Method

Learn how to set up a scan method and acquire data for the sulfa demonstration mix.

3 Qualitative Data Analysis

Learn how to examine chromatograms and spectra to identify sample components. In these exercises, you review data from the sulfa sample you analyzed in Chapter 2, or from a data file that you received with your OpenLab ChemStation software.

4 Set Up and Run a SIM Method

Learn how to set up a selected ion monitoring (SIM) method and acquire data for the sulfa demonstration mix.

5 Set Up and Run a Sequence

Use these exercises to set up an automated sequence for SIM analyses of the sulfa mix at various concentrations, and to acquire data with that sequence.

6 Quantitative Data Analysis

Learn how to analyze data when you need to quantify sample components. These exercises use caffeine data files that you received with your OpenLab ChemStation software.

This page intentionally left blank.

Content

- 1 Prepare for the Analysis 7**
 - Exercise 1. Prepare the LC to run the sample 8**
 - Task 1. Start up ChemStation program 8
 - Task 2. Purge the pump 9
 - Task 3. Prepare the column for the analyses 10
 - Exercise 2. Prepare the samples for the analyses 12**
 - Exercise 3. Check the current MS tune values and adjust if necessary 14**

- 2 Set Up and Run a Scan Method 15**
 - Exercise 1. Set up a full-scan acquisition method 16**
 - Task 1. Enter LC acquisition parameters 16
 - Task 2. Enter MS acquisition parameters 18
 - Task 1. Enter sample information 21
 - Task 2. Acquire the data 22

- 3 Qualitative Data Analysis 25**
 - Exercise 1. Display and manipulate chromatograms 26**
 - Exercise 2. Examine mass spectra 29**
 - Exercise 3. Integrate the chromatogram 33**
 - Exercise 4. Print a report 36**

- 4 Set Up and Run a SIM Method 39**
 - Exercise 1. Set up a SIM acquisition method 40**
 - Task 1. Load the scan method you created previously 40
 - Task 2. Enter MS acquisition parameters 41
 - Exercise 2. Acquire data with the SIM method 43**
 - Task 1. Enter sample information 43
 - Task 2. Acquire the data 45

Content

5	Set Up and Run a Sequence	47
	Exercise 1. Set up a sequence	48
	Task 1. Prepare to create a new sequence	48
	Task 2. Edit sequence parameters	49
	Task 3. Set up the sequence table	51
	Task 4. Set up the sequence output	53
	Exercise 2. Run the sequence	55
6	Quantitative Data Analysis	57
	Exercise 1. Create a method for quantification	58
	Task 1. Create a new method	58
	Task 2. Set up the signal for quantification	59
	Task 3. Integrate the low-level standard	61
	Task 4. Set general calibration parameters	63
	Task 5. Set up the calibration curve	64
	Task 6. Explore options to refine the calibration	68
	Exercise 2. Process a sample and print a report	69

1

Prepare for the Analysis

This chapter presents exercises to help you learn how to:

- Prepare the LC and column for an analysis
- Prepare the samples that you analyze in these exercises

Exercise 1. Prepare the LC to run the sample **8**

Task 1. Start up ChemStation program **8**

Task 2. Purge the pump **9**

Task 3. Prepare the column for the analyses **10**

Exercise 2. Prepare the samples for the analyses **12**

Exercise 3. Check the current MS tune values and adjust if necessary **14**

- Check the tune settings of the MS and adjust if necessary.

Before you start

- Order the sample: Agilent Electrospray LC Demo Sample, *p/n 59987-20033*.
- Order the column: Agilent ZORBAX SB-C18, 2.1 mm x 30 mm, 3.5 μm , *p/n 873700-902*.
 - You may use another similar column, but you may need to adjust the HPLC conditions to obtain good separation.
- Make sure that the electrospray source is installed.
- Read the Quick Start Guide.



Prepare for the Analysis

Exercise 1. Prepare the LC to run the sample

Exercise 1. Prepare the LC to run the sample



For the following tasks, try the steps in the first column. If you need more help, follow the detailed instructions in the center column.

Task 1. Start up ChemStation program

Steps	Detailed Instructions	Comments
1 Open the ChemStation window.	<ul style="list-style-type: none">Click the Control Panel icon on the desktop, and then the ChemStation icon.  	ChemStation is an abbreviation of OpenLab ChemStation.

Task 2. Purge the pump


Use these instructions with the binary and quaternary pumps. See the ChemStation online Help for instructions for the capillary and nanoflow pumps.


Steps	Detailed Instructions	Comments
1 Display the Method and Run Control view.	<ul style="list-style-type: none"> In the view selection area in the lower left, click Method and Run Control. 	
2 Place the pump in standby mode.	<ol style="list-style-type: none"> Click More Binary Pump > Control on the Instrument menu to open the Pump Control dialog box. Select Standby and click OK. 	Alternate method: <ul style="list-style-type: none"> Select Standby from the Pump shortcut menu.
3 Prepare solvents used in these familiarization exercises. <ul style="list-style-type: none"> A – 5 mM ammonium formate in water B – 5 mM ammonium formate in methanol 	<ol style="list-style-type: none"> Into a 1-liter reservoir of HPLC-grade water, add 1 mL of 5 M ammonium formate. Into a 1-liter reservoir of HPLC-grade methanol, add 1 mL of 5 M ammonium formate. 	<ul style="list-style-type: none"> The part number for ammonium formate is G1946-85021. Each ampoule contains 2.2 mL of ammonium formate solution.
4 Replace the solvent bottles with the ones you just prepared.	<ul style="list-style-type: none"> Replace the bottles for channels A and B. 	
5 Open the purge valve.	<ol style="list-style-type: none"> Turn the black purge valve on the front of the pump counter-clockwise two turns. Place the tubing that exits the pump into a 250-mL or larger beaker. 	
6 Enter a flow of 5 mL/min and 50% B, using water in channel A and methanol in channel B.	<ol style="list-style-type: none"> Right-click the pump button. Select Method. Enter the parameters in step 6 and click OK. 	<ul style="list-style-type: none"> Be sure to use HPLC-grade solvents.
7 Turn the pump on and monitor the tubing for bubbles.	<ol style="list-style-type: none"> To turn the pump on, click the green button in the Binary Pump pane. Monitor for bubbles. 	 <ul style="list-style-type: none"> Purge for about 3 minutes to pass 3X the volume for the binary pump. If you wish, you may purge each channel individually first, to ensure that neither is air-locked.

Steps	Detailed Instructions	Comments
8 After the bubbles are gone and the purge is complete, enter a flow of 1 mL/min and 100% B.	<ol style="list-style-type: none"> Right-click the pump button. Select Method. Enter the new parameters in step 8, and click OK. 	
9 Purge a short while longer, and then close the purge valve.	<ol style="list-style-type: none"> Continue to purge for a short while. Close the black valve. 	For more information on purging the pump, see the reference manual that you received with your pump.

Task 3. Prepare the column for the analyses

In the exercises in the next chapters, you analyze a mixture of four sulfonamide compounds. To perform the analyses in the following chapters, you must first condition and equilibrate your column.

Steps	Detailed Instructions	Comments
1 Disconnect the column from the detector and MS.	<ol style="list-style-type: none"> Turn the pump off by clicking the red button in the Binary Pump pane.  Disconnect the column from the detector and MS. Place the open end of the tubing that exits the column into the beaker. 	<ul style="list-style-type: none"> To prevent detector contamination, allow the column effluent to go directly to the waste beaker.
2 Flush the column with 100% methanol at 1 mL/min (5 to 10 min). <ul style="list-style-type: none"> ZORBAX SB-C18, 2.1 mm X 30 mm, 3.5 µm, <i>p/n 873700-902</i> 	<ol style="list-style-type: none"> Turn the pump on. Flow methanol through the column under the conditions used in Task 2, step 8. 	<ul style="list-style-type: none"> The data sheet shipped with the column cartridge recommends that you flush with 20 to 30 column-volumes of 100% methanol (approximately 5 to 7.5 mL).
3 Condition the column as follows, using the solvents made up in Task 2, step 6 : <ul style="list-style-type: none"> Flow rate – 0.4 mL/min 100% B for 1/2 hour 50% B for 1/2 hour 	<ol style="list-style-type: none"> Click Set up Instrument Method on the Instrument menu to open the Setup Method dialog box. Click the Binary Pump tab. Enter the flow rate in step 3. For Solvent B, type 100 and click Apply. Wait 30 minutes. For Solvent B, type 50 and click Apply. Wait 30 minutes. 	<ul style="list-style-type: none"> At a flow rate of 0.4 mL/min, the checkout column should produce about 70 to 80 bar pressure (measured without any fittings at the column exit). If, after you perform these steps, the pump pressure through the column is too high, order a replacement SB-C18 column (<i>p/n 873700-902</i>). If your column is <i>not</i> new, you can reduce the length of time that you condition the column.

Steps	Detailed Instructions	Comments
<p>4 Equilibrate the column at the analysis conditions:</p> <ul style="list-style-type: none"> • 12% B for 1/2 hour at 40 °C 	<p>a For Solvent B, type 12 and click OK. b Click the Column Comp. tab in the Setup Method dialog box. c For Temperature, type 40 and click OK.</p>	<ul style="list-style-type: none"> • While you condition and equilibrate the column, you may complete step 5 in this exercise and then work on the rest of the exercises in this chapter. Be sure to complete step 6 before you go on to the next chapter.
<p>5 While the column equilibrates, set parameters for the MS spray chamber so it can heat and equilibrate as well.</p> <ul style="list-style-type: none"> • Drying gas flow: 8 L/min • Nebulizer pressure: 35 psig • Drying gas temperature: 300 °C • Capillary voltage: 3000 V <p>For Agilent Jet Stream source (LC/MSD XT, 6130B, and 6150B):</p> <ul style="list-style-type: none"> • Sheath Gas Flow: 12 L/min • Sheath Gas Temp: 360°C • Nozzle Voltage: 0 V 	<p>a Right-click the MSD button on the system diagram and select Spray Chamber. </p> <p>b Enter the parameters described in step 5. c Click OK. d Wait 10 minutes before you tune the MS.</p>	
<p>6 Reconnect the column to the DAD and MS.</p>		<ul style="list-style-type: none"> • You can complete “Exercise 3. Check the current MS tune values and adjust if necessary” either with or without the column connected to the DAD and MS, but you <i>do</i> need to reconnect prior to the exercises in Chapter 2, “Set Up and Run a Scan Method.”

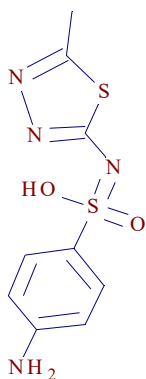
Prepare for the Analysis

Exercise 2. Prepare the samples for the analyses

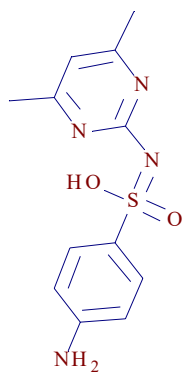
Exercise 2. Prepare the samples for the analyses

In the exercises in the next chapters, you analyze a mixture of four sulfonamide compounds. The Electrospray LC Demo Sample (*p/n* 59987-20033), contains five ampoules with 100 ng/ μ L each of these compounds:

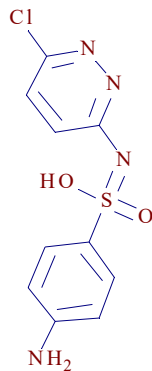
- sulfamethizole (M+H)⁺ = 271
- sulfamethazine (M+H)⁺ = 279
- sulfachloropyridazine (M+H)⁺ = 285
- sulfadimethoxine (M+H)⁺ = 311.



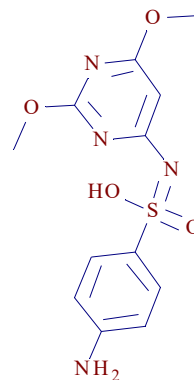
Sulfamethizole



Sulfamethazine



Sulfachloropyridazine




Sulfadimethoxine

To perform the analyses in the following chapters, you must first prepare the sample at various dilutions. The final concentrations will be 1, 5 and 10 ng/ μ L. You will also prepare a solvent blank.

Steps	Detailed Instructions	Comments
<p>1 Prepare a 1:10 dilution of the original sample in a 1-mL autosampler vial.</p> <ul style="list-style-type: none"> Final concentration is 10 ng/μL You will use this sample for the scan analysis in Chapter 2, and for the SIM analyses in Chapter 4 and Chapter 5. 	<p>a Transfer 100 μL of the sulfa mixture into the autosampler vial.</p> <p>b Add 900 μL of 90:10 water:methanol that contains 5 mM ammonium formate (NH₄HCO₂).</p> <p>c Close the vial with a cap.</p>	<ul style="list-style-type: none"> The original sulfa mixture is dissolved in a solvent mixture of 70% water and 30% acetonitrile.
<p>2 Prepare a 1:20 dilution of the original sample in a 1-mL autosampler vial.</p> <ul style="list-style-type: none"> Final concentration is 5 ng/μL You will use this sample for the SIM analysis in Chapter 5. 	<p>a Transfer 50 μL of the sulfa mixture into the autosampler vial.</p> <p>b Add 950 μL of 90:10 water:methanol that contains 5 mM ammonium formate.</p> <p>c Close the vial with a cap.</p>	
<p>3 Prepare a 1:100 dilution of the original sample in a 1-mL autosampler vial.</p> <ul style="list-style-type: none"> Final concentration is 1 ng/μL You will use this sample for the SIM analysis in Chapter 5. 	<p>a Transfer 10 μL of the sulfa mixture into the autosampler vial.</p> <p>b Add 990 μL of 90:10 water:methanol that contains 5 mM ammonium formate.</p> <p>c Close the vial with a cap.</p>	
<p>4 Prepare a solvent blank in a 1-mL autosampler vial.</p> <ul style="list-style-type: none"> You will use this sample for the SIM analysis in Chapter 5. 	<p>a Into the autosampler vial, transfer 990 μL of 90:10 water:methanol that contains 5 mM ammonium formate.</p> <p>b Close the vial with a cap.</p>	

Exercise 3. Check the current MS tune values and adjust if necessary

The MS is very stable and does not need to be tuned very often. You can usually tune just once a month, or once a week at most. You can use the Check Tune program described in this exercise to confirm that the MS is in adjustment.

Steps	Detailed Instructions	Comments
1 Switch to the MSD Tune view.	<ul style="list-style-type: none"> In the view selection area in the lower left, click MSD Tune. 	
2 Select the tune file.	<ol style="list-style-type: none"> In the Select Tune File dialog box, select ATUNES.TUN. Keep the default of Positive Polarity (Standard). Click OK. In the status bar near the top of the MSD Tune view, verify that you see the following: <ul style="list-style-type: none"> Mode is API-ES Source is ESI (electrospray) 	<ul style="list-style-type: none"> Make sure that you use an appropriate calibrant with an appropriate source.
3 Run a Check Tune.	<ul style="list-style-type: none"> In the Tune menu, select Check Tune. Note that Check Tune requires values for comparison that are determined from a previous Autotune. Autotune is normally run during installation. 	<ul style="list-style-type: none"> Check Tune is normally all that you need to do to confirm that the MS settings are correct. If Check Tune indicates a problem with your MS settings, then proceed to step 4 and/or step 5.
4 If Check Tune report suggests that you adjust peak widths or mass axis, then do that.	<ol style="list-style-type: none"> In the Tune menu, select Adjust Mass Peak Width. In the Tune menu, select Calibrate Mass Axis. 	
5 If the Check Tune report shows poor sensitivity, which indicates that your MS settings are significantly out of adjustment, then run a full Autotune.	<ul style="list-style-type: none"> In the Tune menu, select Autotune > Positive Polarity. 	<ul style="list-style-type: none"> The exercises in this manual use only the positive ion mode and standard scan speeds, so it is not necessary to tune for negative polarity or fast scan.

2

Set Up and Run a Scan Method

These exercises show you how to set up a scan data acquisition method for the demonstration sample (sulfa mix) and to acquire data with that method.

Exercise 1. Set up a full-scan acquisition method **16**

Task 1. Enter LC acquisition parameters **16**

Task 2. Enter MS acquisition parameters **18**

Task 1. Enter sample information **21**

Task 2. Acquire the data **22**

The LC parameters that you enter in these exercises are appropriate for the standard Agilent 1100/1200/1260/1290 Series liquid chromatography (LC) systems. You must enter LC parameters that are appropriate for your LC model.

To view the results of these exercises, see **Chapter 3**, “Qualitative Data Analysis.”

Before you start

- Review the *Quick Start Guide*.
- Prepare the LC, column and sample as described in **Chapter 1**, “Prepare for the Analysis.”


For the tasks on the following pages, try the steps on the left without the detailed instructions. If you need more help, follow the detailed instructions on the right.

Exercise 1. Set up a full-scan acquisition method

This exercise changes the default method and saves it as a new method. This exercise consists of the following tasks:


- “**Task 1. Enter LC acquisition parameters**” on page 16
- “**Task 2. Enter MS acquisition parameters**” on page 18

Task 1. Enter LC acquisition parameters

Steps	Detailed Instructions	Comments
1 Display the Method and Run Control view.	<ul style="list-style-type: none"> • In the view selection area in the lower left of the ChemStation window, click Method and Run Control. 	
2 Open the method DEF_LC.M.	<ol style="list-style-type: none"> Select File > Load > Method. If necessary, navigate to C:\Users\Public\Public Documents\Chemstation\1\methods. Select DEF_LC.M and click OK. 	
3 Save the method under a new name, SULFA MS SCAN 1.M.	<ol style="list-style-type: none"> Select File > Save As > Method. In the dialog box, for Name, type SULFA MS SCAN 1.M. Click OK. In the box for Comment for method history, type a comment. Click OK. 	<ul style="list-style-type: none"> • You save the method now with a new name to avoid inadvertently overwriting the default method later.
4 Enter a volume of 1 μ L for the injection.	<ol style="list-style-type: none"> Click Set up Instrument Method in the Instrument menu to open the Setup Method dialog box. Click the HiP Sampler tab. In the Injection volume box, type 1 for a 1-μL injection. 	

Steps	Detailed Instructions	Comments
5 Enter pump parameters.	<p>a Click the Binary Pump tab in the Setup Method dialog box.</p> <p>b Set the parameters as follows: Flow=0.400mL/min StopTime=7.00 min PostTime=3.00min Solvent A=5mM ammonium formate in water Solvent B=5mM ammonium formate in methanol</p>	
6 Set up the gradient timetable: Line 1 Time 1:00, %B=12, Flow=0.4 Line 2 Time 3:00, %B=100, Flow=0.4 Line 3 Time 6:00, %B=100, Flow=0.4 Line 4 Time 7:00, %B=12, Flow=0.4	<p>a Open the Timetable area in the lower part of the tab, click Add, and type the first line.</p> <p>b Click Add and type the second line.</p> <p>c Repeat step b for lines 3 and 4.</p>	
7 Enter a column compartment temperature of 40 °C.	<p>a Click the Column Comp. tab in the Setup Method dialog box.</p> <p>b Select °C.</p> <p>c Type 40.0 for °C.</p>	
8 Enter parameters for the diode-array detector (DAD).	<p>a Click the DAD tab in the Setup Method dialog box.</p> <p>b Enter the parameters shown below:</p> <ul style="list-style-type: none"> • Use Signal A: Wavelength 272nm, Bandwidth 16 nm • Reference Wavelength: 360 nm, Reference Bandwidth 100 nm • Spectrum Store: All in peak • Peakwidth: > 0.10 min <p>c Click OK to close the Setup Method dialog box with the new setpoints.</p>	<ul style="list-style-type: none"> • The DAD is used in this example, but the variable wavelength detector (VWD) may be used analogously.
9 Select Data Acquisition only in the Run Time Checklist.	<p>a Click Run Time Checklist in the Method menu.</p> <p>b Mark the Data Acquisition check box.</p> <p>c Click OK.</p>	<ul style="list-style-type: none"> • While it is common to include Data Analysis in the Run Time Checklist, for these exercises, you will view the results in Chapter 3, "Qualitative Data Analysis."
10 Save the new parameters to the method file, SULFA MS SCAN 1.M.	<p>a Select File > Save > Method.</p> <p>b In the box for Comment for method history, type a comment.</p> <p>c Click OK.</p>	

Task 2. Enter MS acquisition parameters

Steps	Detailed Instructions	Comments
<p>1 Enter parameters for the quadrupole mass spectrometer (MS):</p> <ul style="list-style-type: none"> Signal 1, scan mode, positive polarity Scan range: 100 to 500 Fragmentor: 100 V for the 6120; 125 V for the LC/MSD, LC/MSD XT, 6130 or 6150 Gain: 1.00 Threshold: 150 Stepsize: 0.10 Peakwidth: 0.05 min Scan data storage: Condensed Active signals: 1 only 	<p>a Right-click the MSD button on the system diagram and select Set up MSD Signals.</p> <p>b Enter the parameters described in step 1 and shown in the figure below. Take care to enter the appropriate Fragmentor voltage for your MS model.</p> <p>c Click OK.</p> 	<ul style="list-style-type: none"> To save disk space you usually acquire line spectra (Scan Data Storage = Condensed). However, when you acquire spectra from intact proteins or protein digests/peptides, you must acquire and deconvolute profile spectra. (Scan Data Storage = Full).

Use MSD

StopTime:

FIA Disabled

General

Tune File:

Ion Source:

Peakwidth: min

Cycle Time: sec/cycle

Fast Scan

Time Filter

Scan Data Storage:

Active Signals:

1

2

3

4

Acquisition Parameters

Display EIC Parameters

Signal: Mode:

Polarity: % cycle time:

	Time(min)	On/Off	Mass Range Low	High	Fragmentor	Gain	Threshold	Step size	Speed (u/sec)
1	0.00	<input checked="" type="checkbox"/>	100.00	500.00	125	1.00	150	0.10	867

Set to 100 for 6120
Set to 125 for LC/MSD, LC/MSD XT, 6130, or 6150


Signal: Mode:

Polarity: % cycle time:

	Time(min)	On/Off	Mass Range Low	High	Fragmentor	Gain	Threshold	Step size	Speed (u/sec)
1	0.00	<input checked="" type="checkbox"/>	100.00	1000.00	70	1.00	150	0.10	

Set Up and Run a Scan Method

Task 2. Enter MS acquisition parameters

Steps	Detailed Instructions	Comments
2	<p>Enter parameters for the spray chamber of the ion source:</p> <ul style="list-style-type: none"> Drying gas flow: 9 L/min Nebulizer pressure: 40 psig Drying gas temperature: 300 °C Capillary voltage: 3000 V <p>For Agilent Jet Stream (LC/MSD XT, 6130B, and 6150B):</p> <ul style="list-style-type: none"> Drying gas flow: 7 L/min Nebulizer pressure: 30 psi Drying gas temperature: 350 °C Sheath gas temperature: 360 °C Capillary voltage: 4000 V Nozzle voltage: 2000 V 	<p>a Right-click the MSD button on the system diagram and select Spray Chamber.</p>  <p>b Enter the parameters described in step 2 and shown in the figure below.</p> <p>c Click OK.</p>

MSD Spray Chamber

Method Spray Chamber: API-ES Lamp: ON OFF

Installed Spray Chamber: API-ES

Temperatures, Pressure, and Flow

	Actual	Setpoint	Maximum
Drying Gas Flow (l/min):	9.0	9.0	13.0
Nebulizer Pressure (psig):	40	40	60
Drying Gas Temperature (°C):	313	300	350
Vaporizer Temperature (°C):		N/A	N/A
Sheath Gas Temperature (°C):		N/A	N/A
Sheath Gas Flow (l/min):		N/A	N/A

Parameters

	Positive	Negative
Capillary Voltage (V):	3000	3000
Corona Current (µA):	N/A	N/A
Charging Voltage (V):	N/A	N/A

Time Table

Time (min)	Parameter	Value
------------	-----------	-------

Insert Append Cut Copy Paste

OK Cancel Help

For all models except 6150B with Agilent Jet Stream source

MSD Spray Chamber

Method Spray Chamber: AJS-ES Lamp: ON OFF

Installed Spray Chamber: API-ES

Temperatures, Pressure, and Flow

	Actual	Setpoint	Maximum
Drying Gas Flow (l/min):	9.0	7.0	13.0
Nebulizer Pressure (psig):	40	30	60
Drying Gas Temperature (°C):	352	350	350
Vaporizer Temperature (°C):		N/A	N/A
Sheath Gas Temperature (°C):		150	360
Sheath Gas Flow (l/min):		12.0	12.0

Parameters

	Positive	Negative
Capillary Voltage (V):	4000	1300
Corona Current (µA):	N/A	N/A
Nozzle Voltage (V):	2000	2000


Time Table

Time (min)	Parameter	Value
------------	-----------	-------

Insert Append Cut Copy Paste

OK Cancel Help

For 6150B with Agilent Jet Stream source

Steps	Detailed Instructions	Comments
3 Set up to store the fragmentor voltage throughout the run.	<p>a Right-click the MSD button on the system diagram and select Data Curves.</p> <p>b Select Fragmentor - 1.</p> <p>c Click Add.</p> <p>d Click OK.</p>	
<div style="border: 1px solid gray; padding: 5px;"> <p style="text-align: right; margin: 0;">MSD Data Curves ✕</p> <hr/> <p>Data Curves</p> <div style="display: flex; justify-content: space-between;"> <div style="width: 45%;"> <p>Available</p> <ul style="list-style-type: none"> ChamCur DryingGas <li style="background-color: #e0e0e0;">Fragmentor - 2 Fragmentor - 3 Fragmentor - 4 Gain - 1 Gain - 2 </div> <div style="width: 45%;"> <p>Selected:</p> <div style="border: 1px solid gray; padding: 2px; margin: 5px 0;">Fragmentor - 1</div> <p style="text-align: center; margin: 0;"> <input type="button" value="Add ->"/> <input type="button" value=" < Remove"/> <input type="button" value="Remove All"/> </p> </div> </div> <p style="text-align: right; margin: 10px 0;">Approximate Data Rate: <input type="text" value=".5"/> Hz.</p> <p style="text-align: center; margin: 0;"> <input type="button" value="OK"/> <input type="button" value="Cancel"/> <input type="button" value="Help"/> </p> </div>		
4 Save the method.	<p>a Select Method > Save Method to overwrite the method SULFA MS SCAN 1.M.</p> <p>b In the box for Comment for method history, type a comment.</p> <p>c Click OK.</p>	

Steps	Detailed Instructions	Comments
5 Print the method.	<p>a Click Method > Print Method.</p> <p>b Mark the check boxes as shown in the figure below.</p> <p>c Click Print.</p>	

Print Method: LC_MSD

Select Parts of the Method to be printed: Check All

Miscellaneous

Method Information

Method Audit Trail

Run Time Check List

Instrument/Acquisition

Detectors Injector/AFC Pump/Oven/Chip

Data Analysis

Report Specification Integration Events

Calibration Data LibSearch and Others

Custom Fields


Select Destination for Printout:

Printer File

Now you are ready to acquire data for the sulfa mix with the method you just created. This exercise consists of the following tasks:

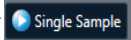

- “**Task 1. Enter sample information**” on page 21
- “**Task 2. Acquire the data**” on page 22

Task 1. Enter sample information

Steps	Detailed Instructions	Comments
1 Display the Single Sample toolbar.	<ul style="list-style-type: none"> • In the top toolbar, click Single Sample. 	

Steps	Detailed Instructions	Comments
2 Display the Sample Information dialog box.	a Click Sample Info in the RunControl menu.	
3 Enter the sample information: <ul style="list-style-type: none"> Operator name Subdirectory: Sulfas Prefix: Sulfa_scan Location: Vial 1 Sample Name: Sulfa 10 ng/μL Comment: Scan familiarization exercise 	<p>a Enter the parameters described in step 3 and shown in the figure below.</p> <p>b Click OK.</p> <ul style="list-style-type: none"> If you select Prefix/Counter, the file names increment automatically from one run to the next. 	

Task 2. Acquire the data

Steps	Detailed Instructions	Comments
1 Place the vial of sulfa sample you prepared at 10 ng/ μ L into position 1 in the autosampler.		<ul style="list-style-type: none"> You prepared this sample in “Exercise 2. Prepare the samples for the analyses” on page 12.
2 Inject the sulfa mix sample.	<ul style="list-style-type: none"> Click Single Sample to start the run. 	<p>This button is present only when you have selected Single Sample mode from the top toolbar.</p> 

Steps	Detailed Instructions	Comments
3 Monitor the total ion chromatogram and the UV chromatogram during data acquisition.	<p>a In the Online Plot window, click Change.</p> <p>b In the list of Available Signals, select DAD A: Signal=272,16 Reference=360,100 and click Add.</p> <p>c In the list of available signals, select MSD: Signal 1 and click Add.</p> <p>d Monitor the MS signal to ensure a stable baseline.</p>	<ul style="list-style-type: none"> If the baseline fluctuation for the MS signal is greater than 10%, the nebulizer and source chamber may require maintenance. See the <i>Agilent InfinityLab LC/MSD Series and 6100 Series LC/MS System Maintenance Guide</i>.
4 Save the signals for the Online Plot window.	<p>a In the Edit Signal Plot dialog box, click Apply to Method.</p> <p>b Save the method.</p>	
5 When the analysis is done, view the results.	<ul style="list-style-type: none"> To view the results, go to the next exercise. 	<ul style="list-style-type: none"> The C18 column may require one or two injections of the sample to be fully conditioned. During these initial injections, everything may be eluted from the column in the void volume. Repeat the process and separation will occur.

This page intentionally left blank.

3

Qualitative Data Analysis

Exercise 1. Display and manipulate chromatograms 26

Exercise 2. Examine mass spectra 29

Exercise 3. Integrate the chromatogram 33

Exercise 4. Print a report 36

This chapter shows you how to analyze data when you need to identify or confirm sample components.

*These exercises use the data file you generated in **Chapter 2**. Alternatively, you can use the sulfa demo data file that you received with the ChemStation software.*



Before you start

- Read the *Quick Start Guide*.
- Set up and run the acquisition method in **Chapter 2**, “Set Up and Run a Scan Method” or that you have the **mssulfas.d** data file in the **MSDEMO** data folder on your system.

For the tasks on the following pages, perform the exercises in the order they appear. Try the steps on the left without the detailed instructions. If you need more help, follow the detailed instructions on the right.

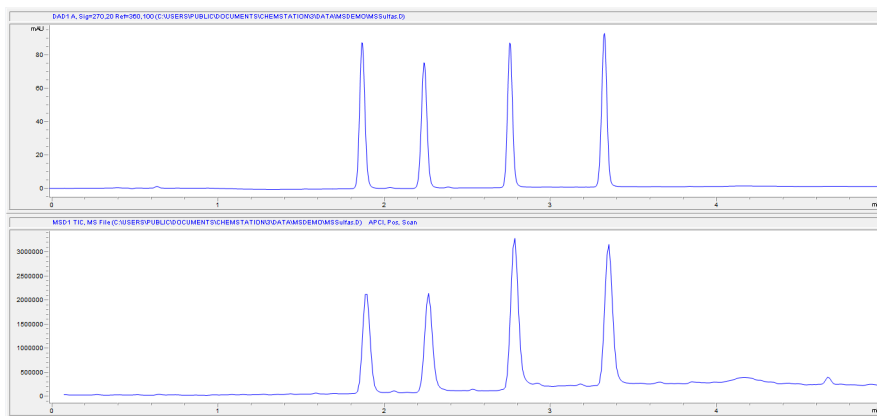
Exercise 1. Display and manipulate chromatograms

In this exercise, you load chromatograms and change the chromatographic display.

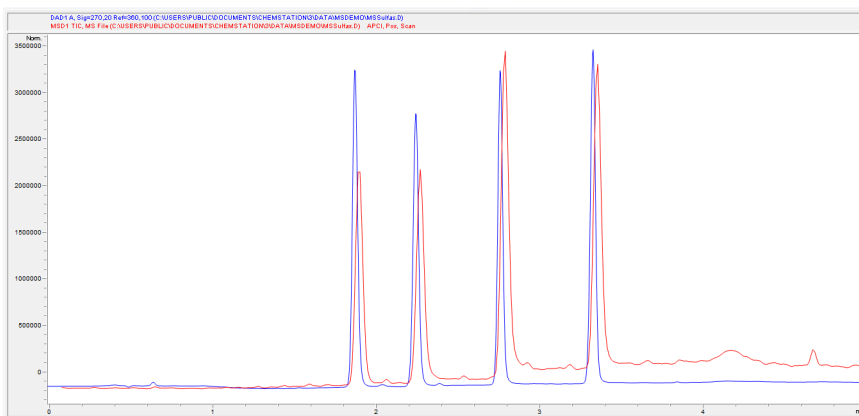
Steps	Detailed Instructions	Comments
1 Display Data Analysis view.	<ul style="list-style-type: none"> In the view selection area of the ChemStation window, click Data Analysis. 	
2 Load the method SULFA MS SCAN 1.M.	<ol style="list-style-type: none"> Select File > Load > Method. Navigate to the folder C:\Users\Public\Public Documents\Chemstation\1\Methods. Select the method file and click OK. 	<ul style="list-style-type: none"> If you just completed the previous exercise, that method is already loaded.
3 Display the Signal Toolset.	<ul style="list-style-type: none"> Click Signal, which is near the center of the window. 	
4 Do one of the following: <ul style="list-style-type: none"> Open the data file, SULFA_SCAN00001.D, which you acquired in Chapter 2. 	<ol style="list-style-type: none"> Select File > Load Signal. Navigate to the appropriate folder, either: <ul style="list-style-type: none"> C:\Users\Public\Public Documents\Chemstation\1\Data\Sulfas, or C:\Users\Public\Public Documents\Chemstation\1\Data\MSDEMO. Select the data file. Set other parameters as shown below and click OK. 	<ul style="list-style-type: none"> If you wish to complete Chapter 4, "Set Up and Run a SIM Method", then you need to process the data file you generated in Chapter 2. You need the report from that data file to set up your SIM groups.

Steps	Detailed Instructions	Comments
<p>File name: MSSulfas.D</p> <p>Ca1Cal01.D Ca1Cal02.D Ca1Cal03.D Ca1Cal04.D Ca1Cal05.D FIA.D LoadTest.D MS3frag.D MSAlign.D MSPosNeg.D MSPurity.D MSScnSM.D MSSulfas.D Prot_FIA.D Protein.D</p> <p>File Information...</p> <p><input type="checkbox"/> Load using Signal Details</p> <p>Signal Information Spectra: dad1: 336 Spectra MSD: 427 Cond.</p> <p><input type="checkbox"/> Integrate after load <input checked="" type="checkbox"/> Load from BSB <input type="checkbox"/> Integrate and print report after load</p>	<p>Folders: c:\...</p> <p>Public Documents ChemStation 3 Data MSDemo</p> <p>Drives: c: Windows</p> <p>Signal Details...</p> <p>Signals: DAD1 A, Sig=270.20 Ref=360.100 MSD1 TIC, MS File, Pos, Scan</p>	<p>OK</p> <p>Cancel</p> <p>Help</p> <p>Short <<</p> <p>Inst. Curves...</p> <p>Network...</p>

- 5 Verify that you see the DAD and MS chromatograms.
- Check that you see a display that is similar to the one shown below.
 - Verify that you see the DAD signal in the top chromatogram.
 - Confirm that you see the MSD signal in the bottom chromatogram.



Steps	Detailed Instructions	Comments
6	<p>Change the chromatogram view so that the MS and UV chromatograms are overlaid in the display.</p> <p>a In the Signal Toolset near the center of the window, click Display Overlaid Signals.</p> <p>b Check that you see the overlaid chromatograms, as shown below.</p> <p>c Click Separate Signals.</p>	<p>The button in step a is also available in the Graphics Toolset, but in that toolset it toggles overlaid/separate. You click the button shown above to turn on the display of the Graphics Toolset.</p>



7	<p>Remove the DAD signal from the display.</p> <p>a In the Navigation Table, click + to display more information.</p> <p>b Under the Signals tab, double-click the signal labeled MSD1 TIC.</p> <p>c When you see the message about the method, click OK.</p> <p>d Verify that the DAD window is gone and only the TIC is displayed.</p>	<p>If you do not see the Navigation Table shown below, in the top toolbar, click the button shown above.</p>
---	--	--

Overlay	Type	Date Time	Operator	Vial	Reference	Data File	Sample Name	Acq. Method
<input type="checkbox"/>		10/29/1997 11:11:11...	C. Miller	1		MSSulfas.D	sulfa drug mix	MSPURITY.M


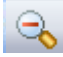
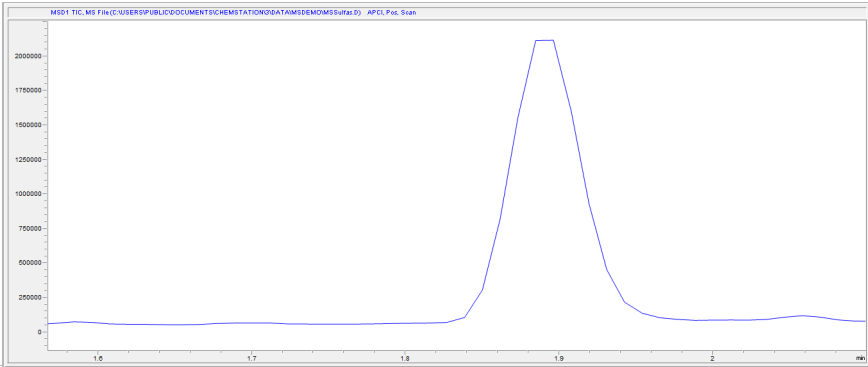
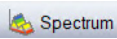
a




Overlay	Type	Date Time	Operator	Vial	Reference	Data File	Sample Name	Acq. Method
<input type="checkbox"/>		10/29/1997 11:11:11...	C. Miller	1		MSSulfas.D	sulfa drug mix	MSPURITY.M
Signals General Info Instrument Curves								
Description		Load?						
DAD 1 A, Sig=270,20 Ref=360,100		<input checked="" type="checkbox"/>						
MSD1 TIC, MS File, Pos, Scan		<input checked="" type="checkbox"/>						

b

Exercise 2. Examine mass spectra

In this exercise, you learn to display mass spectra. You choose background (reference) spectra that you can later subtract from the spectra of the peaks of interest. You learn how to display a single spectrum and an averaged spectrum for a peak.

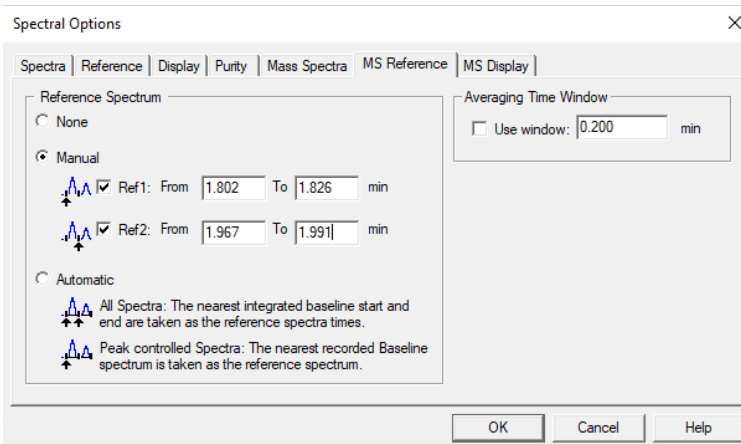
Steps	Detailed Instructions	Comments
1 Zoom in on the first peak in the chromatogram.	<p>a Click Zoom In.</p> <p>b Use the mouse pointer to draw a rectangle around the peak. Take care to include the chromatographic baseline.</p> <p>c Check that your peak looks similar to the one below.</p> <p>d Note the width of the peak at half height. You will need this information to set up the SIM analysis in Chapter 4.</p>	<p>• If you want to try again, you can zoom back out. Do one of the following:</p> <ul style="list-style-type: none"> • Double-click the chromatogram window. • Click Zoom Out.
		 
		
2 Display the Spectrum Toolset.	<p>a Click Spectrum, which is near the center of the window.</p> <p>b If there is not room under your chromatogram window to display spectra, use your mouse pointer to reduce the height of the chromatogram window.</p>	

Steps	Detailed Instructions	Comments
3 Get the first reference spectrum, to the left of the peak.	<p>a To select the first reference spectrum, click the button that is highlighted here.</p>  <p>b In the chromatogram window, do one of the following at the chromatographic baseline just before the peak:</p> <ul style="list-style-type: none"> • Click to select a single spectrum. • Click and drag to select an average spectrum. 	
4 Get a second reference spectrum, to the right of the peak.	<p>a To select the second reference spectrum, click the button that is highlighted here.</p>  <p>b In the chromatogram window, do one of the following at the chromatographic baseline just after the peak:</p> <ul style="list-style-type: none"> • Click to select a single spectrum. • Click and drag to select an average spectrum. 	
5 View your reference spectra.	<p>a If you cannot see the spectra, adjust the size and location of the window labeled Reference Mass Spectrum(a).</p> <p>b Note the two background spectra — one before the peak and one after.</p>	
6 Set the spectral options to do manual background subtraction.	<p>a Click the button to display the Spectral Options dialog box.</p>  <p>b Click the MS Reference tab.</p> <p>c Under Reference Spectrum, click Manual.</p> <p>d Mark the check boxes for Ref1 and Ref2. Note that the time ranges of the reference spectra that you just selected are specified there.</p> <p>e Click OK.</p>	<ul style="list-style-type: none"> • The spectral options apply to all subsequent spectra until you change the options. • If the chromatographic baseline changes over the course of the run, select new reference spectra that are close in time to each peak of interest. • Near the center of the Data Analysis window, you can view and change the setting for background subtraction. <div style="border: 1px solid gray; padding: 2px; width: fit-content; margin-top: 10px;">MSD, Manual Reference ▾</div>


Steps

Detailed Instructions

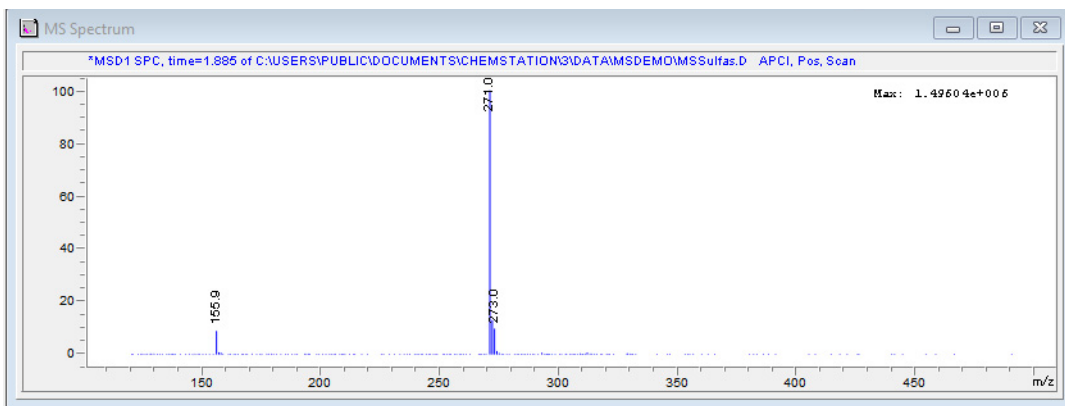
Comments



7 Get a single background-subtracted spectrum for the first LC peak.


- Click the button to get a mass spectrum at any time point. 
- In the chromatogram window, click somewhere on the peak to get the spectrum.
- If necessary for easier viewing, adjust the size and location of the window labeled **MS Spectrum**.
- Verify that the spectrum is similar to the one shown below.

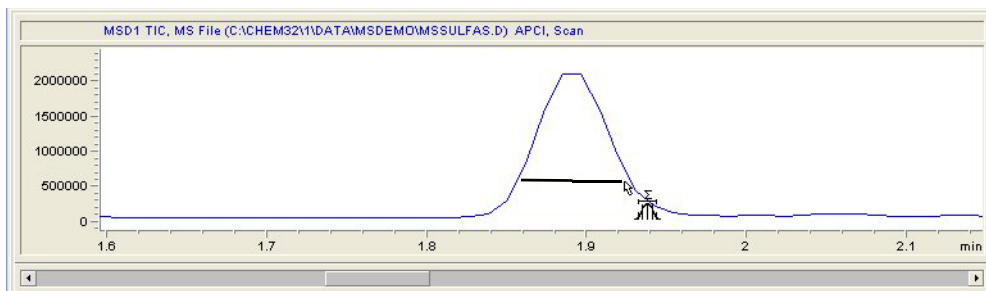
- Under the conditions used to acquire the demo data file (mssulfas.d), the compounds elute in the following order:
Sulfamethizole, $m/z = 271$
Sulfachloropyridazine, $m/z = 285$
Sulfamethazine, $m/z = 279$
Sulfadimethoxine, $m/z = 311$
- Depending on the organic mobile phase and the modifiers, the elution order for the 279 and 285 may change.



Qualitative Data Analysis

Exercise 2. Examine mass spectra


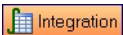
Steps	Detailed Instructions	Comments
8 Get a average background-subtracted spectrum for the first LC peak.	<p>a Click the button to get an averaged mass spectrum.</p> <p>b In the chromatogram window, click and drag the mouse across the peak, as shown below.</p> <p>c View the average spectrum in the window labeled MS Spectrum.</p>	 <ul style="list-style-type: none">When a chromatographic peak consists of a single compound, an average spectrum is usually more accurate.

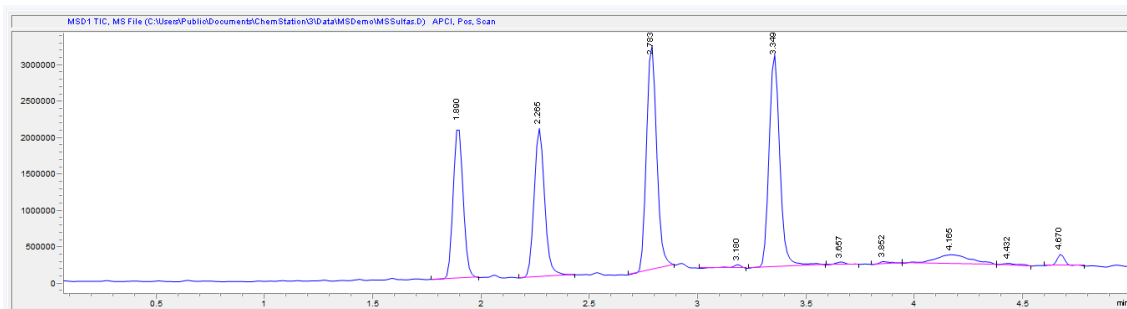


- 9 Be sure to see step 6 in "Exercise 3. Integrate the chromatogram" for an easier, faster way to display spectra.

Exercise 3. Integrate the chromatogram

In this exercise, you learn to set integration events and integrate the chromatogram. Even if you do not care about quantitation, integration helps locate peaks for other purposes. For example, after integration, mass spectra of each peak can be printed with a report.

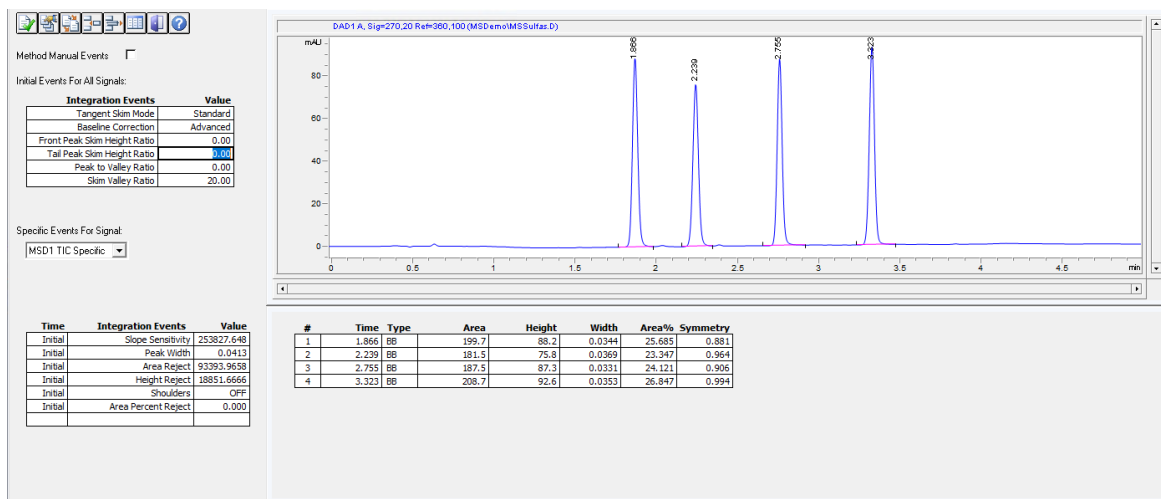
Steps	Detailed Instructions	Comments
1 Display the total ion chromatogram in its entirety.	<p>a Minimize any spectral windows that hide the chromatogram window.</p> <p>b Click Zoom Out.</p>	
2 Display the Integration Toolset.	<p>• Click Integration, which is near the center of the window.</p>	
3 Integrate the chromatogram.	<p>a Click Auto Integrate, which is near the center of the window.</p> <p>b Verify that the results are similar to those shown below.</p>	<p>• Auto Integrate estimates initial integration parameters.</p> <p>• If you do not see the retention times, in the graphics tools, click the button to display retention times.</p> <p>• If you do not see the pink integration baseline, in the graphics tools, click the button to display baselines.</p>



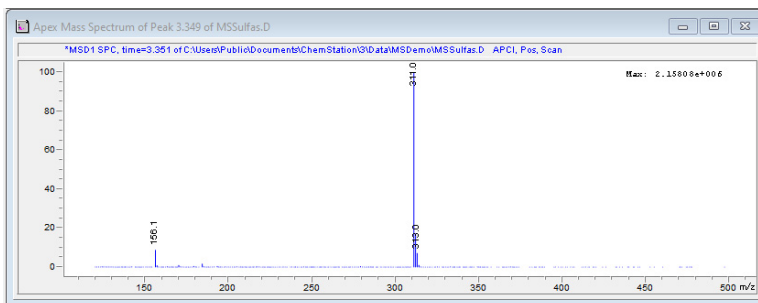
Qualitative Data Analysis

Exercise 3. Integrate the chromatogram

Steps	Detailed Instructions	Comments
4	<p>Adjust the integration parameters to get only four integrated peaks.</p> <p>a Click Edit/Set Integration Events Table.</p> <p>b In the Integration Events table, for Baseline Correction, select Advanced.</p> <p>c For Height Reject, type 500000.</p> <p>d Click Integrate current Chromatogram.</p> <p>e Verify that your results are similar to those shown below.</p>	<p>For detailed information about integration events, see <i>OpenLab ChemStation Concepts and Workflows</i>.</p>



Steps	Detailed Instructions	Comments
5 Save the integration events to the method in memory.	<ul style="list-style-type: none"> Click the button to exit and save the integration results. 	<ul style="list-style-type: none"> To save the events to the method on disk, you also need to save the method to disk, as described in step 3 on page 37.
6 Use the integrated chromatogram as the basis for a faster way to display background-subtracted spectra.	<ul style="list-style-type: none"> a Click Spectrum. b Click the buttons to display the Spectral Options dialog box. c Click the MS Reference tab. d Under Reference Spectrum, click Automatic. e Click OK. f Click the button to get a mass spectrum at the peak apex. g In the chromatogram window, click somewhere on the <i>fourth</i> peak to get the spectrum. h Verify that the spectrum is similar to the one shown below. 	<ul style="list-style-type: none"> When you set Reference Spectrum to Automatic, the software automatically takes the reference spectra for each peak, as described in the Spectral Options dialog box. The button to get the mass spectrum at the peak apex is available only if you have integrated the chromatogram. No matter where you click on the peak, it gets the spectrum at the apex. With this tool, you may not need to zoom in on the chromatogram to achieve a precise location for the spectrum.



Exercise 4. Print a report

In this exercise, you print a report, which you will use in **Chapter 4**, “Set Up and Run a SIM Method.”

Steps	Detailed Instructions	Comments
1	Specify the LCMS Qualitative report style, with the report printed to the screen.	<p>a Select Report > Specify Report.</p> <p>b In the Specify Report dialog box, under Destination, mark the check box for Screen.</p> <p>c For Report Style, select LCMS Qualitative.</p> <p>d Check that other settings are as shown below.</p> <p>e Click OK.</p>

Specify Report: LC_MSD ×

Reporting settings Quantitation settings

Use Intelligent Reporting Use Classic Reporting

Style

Report style: LCMS Qualitative Report Style Options...

Quantitative result sorted by: Signal

Sample info on each page Add fraction table and ticks

Add chromatogram output Add summed peaks table

Add sample custom fields to sample info Add compound custom fields

Report layout for uncalibrated peaks

Separately With calibrated peaks Do not report

Chromatogram output

Portrait

Landscape

Multi-page

1 Pages

Signal options...

Size (% of page)

Time: 100

Response: 40

Destination

Printer

Screen

File

File setting

File prefix: Report

PDF CSV HTM DIF

Unique PDF file name TXT XLS EMF

OK Cancel Help

Steps	Detailed Instructions	Comments
2 Print the report.	<ul style="list-style-type: none"> a Select Report > Print Report. b After a short wait, view the Report window. c Verify that page 1 of the report contains header information and an integrated chromatogram. d At the bottom of the report window, click Next. e Verify that page 2 of the report shows extracted ion chromatograms and a mass spectrum of the first chromatographic peak. f Continue to click Next to view results for the three additional chromatographic peaks. g At the bottom of the report window, click Print. This prints a hard copy of the report. h At the bottom of the report window, click Close. 	<ul style="list-style-type: none"> • If you wish to complete Chapter 4, "Set Up and Run a SIM Method", then save the hard copy and refer to it when you set up your SIM groups. • The extracted ion chromatograms are indicators of peak purity; if the retention times fail to coincide, the peak likely represents more than one compound.
3 Save the method.	<ul style="list-style-type: none"> a Select File > Save > Method to overwrite the method SULFA MS SCAN 1.M. b In the box for Comment for method history, type a comment. c Click OK. 	<ul style="list-style-type: none"> • You save the method now so that your integration parameters, spectral display options, report settings, and other data analysis settings become part of the method.

This page intentionally left blank.

4

Set Up and Run a SIM Method

Exercise 1. Set up a SIM acquisition method 40

Task 1. Load the scan method you created previously 40

Task 2. Enter MS acquisition parameters 41

Exercise 2. Acquire data with the SIM method 43

Task 1. Enter sample information 43

Task 2. Acquire the data 45

These exercises show you how to set up a data acquisition method that uses selected ion monitoring (SIM). You set up the method for the demonstration sample (sulfa mix) and then run the sample with that method.

To set up the SIM method, you modify the scan method that you created in **Chapter 2**. To set up the SIM acquisition, you need the following for each of the four sulfa compounds:

- The LC retention time
- The masses of ions in the spectrum

You get that information from the report you generated in **Chapter 3**.

Before you start


- Complete the previous exercises in this manual.

Exercise 1. Set up a SIM acquisition method


In this exercise, you start with your existing scan method and modify it for SIM analysis. You keep the same LC conditions and modify only the MS conditions. This exercise consists of the following tasks:

- “**Task 1. Load the scan method you created previously**” (below)
- “**Task 2. Enter MS acquisition parameters**” on page 41

Task 1. Load the scan method you created previously

Steps	Detailed Instructions	Comments
1 Display Method and Run Control view.	<ul style="list-style-type: none"> • In the view selection area of the OpenLab ChemStation window, click Method and Run Control. 	
2 Open the method SULFA MS SCAN 1.M.	<ol style="list-style-type: none"> Select File > Load > Method. If necessary, navigate to C:\Users\Public\Public Documents\Chemstation\1\Methods. Select SULFA MS SCAN 1.M and click OK. 	
3 Save the method under a new name, SULFA MS SIM 1.M.	<ol style="list-style-type: none"> Select File > Save As > Method. In the dialog box, for Name, type SULFA MS SIM 1.M. Click OK. In the box for Comment for method history, type a comment. Click OK. 	<ul style="list-style-type: none"> • You save the method now to avoid inadvertent overwrites of your scan method.

Task 2. Enter MS acquisition parameters

Steps	Detailed Instructions	Comments
<p>1 Enter the chromatographic peak width for the SIM analysis.</p>	<p>a Right-click the MSD button on the system diagram and select Set up MSD Signals.</p>  <p>b When the Set Up MSD Signals dialog box is displayed, type 0.05 for Peakwidth.</p>	<ul style="list-style-type: none"> The peak width is an important setting. It is used to calculate the appropriate SIM dwell times to deliver sufficient points across a chromatographic peak to give good quantitation. Peak width is defined as the full width at half maximum (FWHM), the width at 50% of the peak height.
<p>2 Set up the first SIM ions using the masses (to the nearest 0.1) that you observed in the spectra from your scan analysis:</p> <ul style="list-style-type: none"> Sulfamethizole: Time 0, SIM Ions 271 and 156. 	<p>a Under MSD Signal Settings, Signal 1, for Mode, select SIM.</p> <p>b In the table, for Fragmentor, type 200.</p> <p>c In the table, change Group 1 to Sulfamethizole, and for SIM Ion, refer to the spectrum on your printout and type the mass (to the nearest 0.1) for the 271 ion.</p> <p>d Click Add Ion, and type the mass for the sulfamethizole 156 ion.</p>	<ul style="list-style-type: none"> In this example, each SIM group includes a pseudo-molecular ion and one fragment ion for confirmation. Note that the figure below does not show the fourth sulfa drug.
<p>3 Set up the remaining SIM ions, using the masses (to the nearest 0.1) that you observed in the spectra from your scan analysis:</p> <ul style="list-style-type: none"> Sulfachloropyridazine: Time 1.3, SIM Ions 285, 287, and 156. Sulfamethazine: Time 2.3, SIM Ions 279 and 186. Sulfadimethoxine: Time 3.3, SIM Ions 311 and 156. 	<p>a Click Add Grp, and type the name, start time and mass (approximately 285) for sulfachloropyridazine.</p> <p>b Click Add Ion, and type the mass for the sulfachloropyridazine 156 ion.</p> <p>c Click Add Ion, and type the mass for the sulfachloropyridazine 287 ion.</p> <p>d Repeat these steps until you have entered two or three ions for each of the remaining compounds.</p> <p>e Click OK to close the Set Up MSD Signals dialog box.</p>	<ul style="list-style-type: none"> Alternatively, instead of making separate groups for each compound as described here, <i>all</i> of the SIM ions could be entered into "Group 1", which could be re-named "Sulfonamides". The first SIM group can contain up to 100 ions. You may need to adjust the start time for each SIM group. Refer to your printout from Chapter 3 to determine a start time so that each group change occurs about midway between the chromatographic peaks. If the retention time difference between sulfachloropyridazine and sulfamethazine is less than 0.3 minutes, merge these ions into one group. The sulfachloropyridazine additionally includes the chlorine isotope at m/z 287.

Steps

Detailed Instructions

Comments

Set Up MSD Signals

MSD Control
 Use MSD
 StopTime: noLimit
 FIA Disabled
 Tune File: skunes.tun
 Ion Source: API-ES
 Peakwidth: 0.050 min
 Cycle Time: 0.30 sec/cycle
 Fast Scan
 Time Filter
 Scan Data Storage: Condensed
 Active Signals:
 1 SIM_Pos
 2
 3
 4
 Acquisition Parameters
 Display EIC Parameters
 Advanced...

MSD Signal Settings

Signal: 1 Mode: SIM SIM on Sample Target Masses
 Polarity: Positive % cycle time: 100.0

Time(min)	On/Off	Group	SIM Ion	Frag-mentor	Gain	Dwell (msec)	%Rel Dwell	Compound Name	ISTD
1	0.00	✓	Sulfamethiz	156.10	150	1.00	144	50.0	
1				271.00	150		144	50.0	
2	1.00	✓	Sulfamethiz	156.10	150	1.00	95	33.3	
2				285.00	150		95	33.3	
2				287.00	150		95	33.3	
3	2.00	✓	Sulfamethiz	186.10	150	1.00	144	50.0	

Sort Add Ion Add Grip Cut Copy Paste

Signal: 2 Mode: Scan
 Polarity: Positive % cycle time:

Time(min)	On/Off	Mass Range Low High	Frag-mentor	Gain	Thres-hold	Step size	Speed (u/sec)	
1	0.00	✓	100.00	1000.00	70	1.00	150	0.10

Sort Inset Append Cut Copy Paste

OK Cancel Help

4 Save the method.


- a Select **Method > Save Method** to overwrite the method **SULFA MS SIM 1.M**.
- b In the box for **Comment for method history**, type a comment.
- c Click **OK**.

Exercise 2. Acquire data with the SIM method

Now you are ready to acquire data for the sulfa mix with the method you just created. This exercise consists of the following tasks:

- “**Task 1. Enter sample information**” on page 43
- “**Task 2. Acquire the data**” on page 45

Task 1. Enter sample information

Steps	Detailed Instructions	Comments
1 Display the Single Sample toolbar.	<ul style="list-style-type: none">• In the top toolbar, click Single Sample.	
2 Display the Sample Information dialog box.	<ul style="list-style-type: none">a Click Sample Info in the RunControl menu.	

Set Up and Run a SIM Method

Task 1. Enter sample information

Steps	Detailed Instructions	Comments
<p>3 Enter the sample information:</p> <ul style="list-style-type: none"> Subdirectory: Sulfas Prefix: Sulfa_SIM Location: Vial 1 Sample Name: Sulfas 10 ng/μL Comment: SIM familiarization exercise 	<p>a Enter the parameters described in step 3 and shown in the figure below.</p> <p>b Click OK.</p>	<ul style="list-style-type: none"> If you select Prefix/Counter, the file names increment automatically from one run to the next.

Sample Info: LC_MSD

Operator name: SYSTEM

Data file

Path: C:\Users\Public\Documents\ChemStation\3\Data\ Subdirectory: Sulfas

Name Pattern

Signal 1: Sulfa_SIM:Counter: Sulfas_SIM000001.D Counter: 1 Length: 6 Increment: 1

Injection Source: as Method

Error Method

Sample parameters

Sample Location: 1 (blank run if no entry) Plate ID:

Sample name: Sulfas 10 ng/ μ L Sample amount: 0 Multiplier: 1 Dilution: 1

Run Type: Sample

Calibration Settings

Level: 1 Update RF: No Update Update RT: No Update

Lims IDs

1: 2: 3:

ISTD Amounts

1: 0 2: 0

3: 0 4: 0

5: 0 6: 0

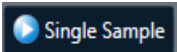

7: 0 8: 0

Comment: Scan familiarization exercise

Target masses:

Custom Fields ... Number of runs: 1 Run Method OK Cancel Help

Task 2. Acquire the data

Steps	Detailed Instructions	Comments
1 Place the vial of sulfa sample you prepared at 10 ng/μL into position 1 in the autosampler.		<ul style="list-style-type: none"> You prepared this sample in “Exercise 2. Prepare the samples for the analyses” on page 12.
2 Inject the sulfa mix sample.	<ul style="list-style-type: none"> Click Single Sample. 	<p>This button is present only when you have selected Single Sample mode from the top toolbar. </p>
3 Monitor the total ion chromatogram and the UV chromatogram during data acquisition.	<ul style="list-style-type: none"> a Activate the Online Plot window. b Monitor the MS signal to ensure a stable baseline. 	<ul style="list-style-type: none"> If the baseline fluctuation for the MS signal is greater than 10%, the nebulizer and source chamber may require maintenance. See the <i>Agilent InfinityLab LC/MSD Series and 6100 Series LC/MS Maintenance Guide</i>.
4 When the analysis is done, view the results.	<ul style="list-style-type: none"> a Display Data Analysis view. b Load the data file you just created. c Examine the DAD and MS chromatograms. 	<ul style="list-style-type: none"> If you need help, follow the general procedure in “Exercise 1. Display and manipulate chromatograms” on page 26 in Chapter 3.

This page intentionally left blank.

5

Set Up and Run a Sequence

Exercise 1. Set up a sequence 48

Task 1. Prepare to create a new sequence 48

Task 2. Edit sequence parameters 49

Task 3. Set up the sequence table 51

Task 4. Set up the sequence output 53

Exercise 2. Run the sequence 55

These exercises show you how to set up a sequence for the SIM analysis of the demonstration sample (sulfa mix), and to acquire data with that sequence.

In the sequence, you run the sulfa mix at three concentrations: 1, 5 and 10 ng/μL. You also run a solvent blank.

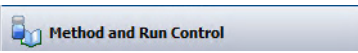

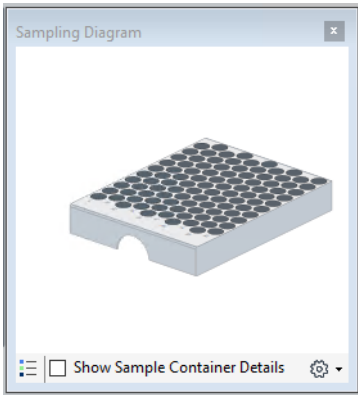
Before you start

- Read the *Quick Start Guide*.
- Complete the previous exercises in this manual.

For details about sequences, see the automation chapter in *OpenLab ChemStation: Concepts and Workflows*.

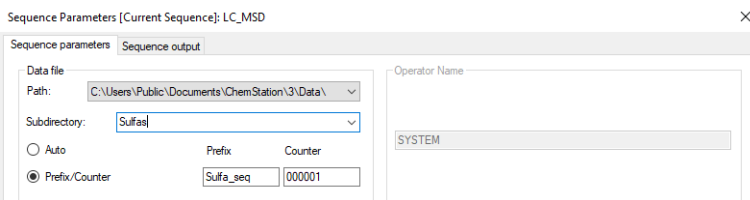
Exercise 1. Set up a sequence

Task 1. Prepare to create a new sequence

Steps	Detailed Instructions	Comments
1 Display Method and Run Control view.	<ul style="list-style-type: none"> In the view selection area of the OpenLab ChemStation window, click Method and Run Control. 	
		
2 Display the Sequence Toolset.	<ul style="list-style-type: none"> In the top toolbar, click the button to display the Sequence Toolset. 	
3 Display the Autosampler Tray diagram.	<ul style="list-style-type: none"> Click Sampling Diagram in the View menu. 	
		
4 Initiate setup of a new sequence.	<ul style="list-style-type: none"> Select Sequence > Sequence Table. 	
5 Save the sequence under a new name, SULFA MS SIM 1.S	<ol style="list-style-type: none"> Select Sequence > Save Sequence Template As. For Name, type SULFA MS SIM 1.S. Click OK. 	

Task 2. Edit sequence parameters

Steps	Detailed Instructions	Comments
1 Open Sequence Parameters dialog box.	<ul style="list-style-type: none"> Click Sequence > Sequence Parameters. 	<ul style="list-style-type: none"> The sequence parameters are settings that are common to all the samples in the sequence.
2 Enter the sequence parameters for Data File.	<p>a Enter the following parameters, shown in step 1.</p> <ul style="list-style-type: none"> Subdirectory: Sulfas Prefix: Sulfa_seq 	<ul style="list-style-type: none"> To avoid overwrite of data files, type a new subdirectory for each sequence. The directory will be created if it doesn't already exist on your computer. Unique file names are automatically created for each data file within the subdirectory.



Steps	Detailed Instructions	Comments
3 Enter the rest of the sequence parameters:	<p>a Enter the following parameters shown in the figure below.</p> <ul style="list-style-type: none"> Parts of methods to run: According to Runtime Checklist Wait: 10 minutes after loading a new method Shutdown: STANDBY Not Ready Timeout: 15 minutes Sequence Comment: Sequence familiarization exercise <p>b Click OK.</p>	<ul style="list-style-type: none"> If you wanted to run only reprocessing (data analysis), you would set that in Part of methods to run. The Wait allows the instrument to equilibrate when a new method is loaded.

Post-Sequence Command/Macros are a convenient way to turn off lamps, pumps, etc. The command or macro is run at the end of the sequence or in the event of an error.

Two examples of Post-Sequence Command/Macros are:

- MSSetState is a command that can change the MS state to standby. See the online Help for commands.
- SHUTDOWN.MAC is a macro that will shut down the system, but you must customize it before using it.

Task 3. Set up the sequence table

Steps	Detailed Instructions	Comments
<p>1 Set up the sequence table to:</p> <ul style="list-style-type: none"> Run duplicate injections of a blank. Run duplicate injections of the sulfa mix at three concentrations: 1, 5 and 10 ng/μL. Use the method SULFA MS SIM 1.M, that you created in Chapter 4, "Set Up and Run a SIM Method". 	<p>a Click Sequence > Sequence Table.</p> <p>b In the first line, type these values:</p> <ul style="list-style-type: none"> Sample Location = 1 Sample Name = ng/ul sulfas Method Name = SULFA MS SIM 1 Inj/Loc = 2 Sample Type = Sample <p>c Select Line 1.</p> <p>d Click Insert/Filldown Wizard.</p> <p>e Complete the Insert/Filldown Wizard as shown in the next figure.</p>	<ul style="list-style-type: none"> In this step, you set up the parts of the sequence table that are common to all the samples. You will specify the sample names later in this exercise. There are a number of ways to add samples to the sequence table. This exercise illustrates just one of the ways – use of the Insert/Filldown Wizard.

Filldown Options: LC_MSD

Define rules to be used for filldown.

Sample Locations

Increment starting from: 2 by: 1

Rectangular increment starting from: 1 to:

Increment location only for sample type: "Sample"

Use 1

No filldown

Increment pattern

Sample Name

Detect counter and increment

Append counter and increment starting at: 0001

Increment only for sample type: "Sample"

Use:

No filldown

Same name for identical sample location

OK Cancel Help

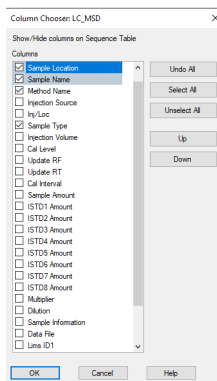
Set Up and Run a Sequence

Task 3. Set up the sequence table

Steps	Detailed Instructions	Comments
2	<p>View the sequence table that you have created so far.</p> <p>a Compare your table with the one below.</p> <p>b Note any differences, such as columns that are included and column widths.</p>	<ul style="list-style-type: none"> Your results will likely differ, but in the next step you can recreate the table format below.

Line	Sample Location	Sample Name	Method Name	Inj/Loc	Sample Type
1	1	Blank	SULFA MS SIM 1	2	Sample
2	2	1 ng/ul sulfas	SULFA MS SIM 1	2	Sample
3	3	5 ng/ul sulfas	SULFA MS SIM 1	2	Sample
4	4	10 ng/ul sulfas	SULFA MS SIM 1	2	Sample

3	<p>(Optional) Customize the sequence table to match the format in step 2.</p> <p>a Click the Column Chooser button.</p> <p>b Click Unselect All to clear all check boxes.</p> <p>c Click OK.</p>	 <ul style="list-style-type: none"> For descriptions of any columns you removed, see the online Help.
---	---	---



4	<p>Type the following sample names into the table:</p> <ul style="list-style-type: none"> Vial 1 – blank Remaining vials – sulfa mix at 1, 5 and 10 ng/μL <p>a Change the Sample Name for each sample, as shown below.</p> <p>b Click OK.</p>
---	---

Line	Sample Location	Sample Name	Method Name	Inj/Loc	Sample Type
1	1	Blank	SULFA MS SIM 1	2	Sample
2	2	1 ng/ul sulfas	SULFA MS SIM 1	2	Sample
3	3	5 ng/ul sulfas	SULFA MS SIM 1	2	Sample
4	4	10 ng/ul sulfas	SULFA MS SIM 1	2	Sample

5	<p>Save the sequence.</p> <p>a Click Save Sequence in the Sequence toolset.</p>	
---	--	---


Task 4. Set up the sequence output

Steps	Detailed Instructions	Comments
1 Set up the sequence to print a short sequence summary to the printer.	<p>a Click Sequence > Sequence Parameters > Sequence Output.</p> <p>b Mark the check box for Print sequence summary report.</p> <p>c Mark the check box for Report to printer.</p> <p>d Click Setup.</p> <p>e Fill in the dialog box as shown below.</p> <p>f Click OK in the Sequence Summary Parameters dialog box.</p> <p>g Click OK in the Sequence Parameters dialog box.</p>	<ul style="list-style-type: none"> In addition to the sequence summary report, you can print individual sample reports, as specified in your method. (You do not print individual reports in this exercise.) For details about sequence reports, see the chapter on ChemStation reports in <i>OpenLab ChemStation: Concepts and Workflows</i>. The setup shown in the dialog box below prints the simplest summary report.

Sequence Summary Parameters [Current Sequence]: LC_MSD

The screenshot shows the 'Sequence Summary Parameters' dialog box with the 'Reports' tab active. Under 'Activate Report', there is a list of reports with checkboxes. Report 9, 'Summary', is checked. To the right, under 'Style', there are three dropdown menus. The first two are set to 'Standard Statistics' and the third, corresponding to report 9, is set to 'Sample Summary'.

Activate Report	Style:
<input type="checkbox"/> 1. One Page Header	
<input type="checkbox"/> 2. Configuration	
<input type="checkbox"/> 3. Sequence	
<input type="checkbox"/> 4. Logbook	
<input type="checkbox"/> 5. Methods	
<input type="checkbox"/> 6. Analysis Reports	
<input type="checkbox"/> 7. Statistics calib. runs	Standard Statistics
<input type="checkbox"/> 8. Statistics sample runs	Standard Statistics
<input checked="" type="checkbox"/> 9. Summary	Sample Summary

Steps	Detailed Instructions	Comments
2 Save the sequence.	<ul style="list-style-type: none"> Click Save Sequence in the Sequence toolset. 	
3 Print the sequence.	<ol style="list-style-type: none"> Select Sequence > Print Sequence Template. Mark the check boxes as shown in the figure below. Click Print. 	<p>If you click the Print All button, you print all the parts of the sequence rather than the items you just specified.</p>

Print Sequence: LC_MSD

Select Parts of the Sequence to be printed:

Sequence Parameters
 Chemstore Transfer Settings

Sequence Table

Sample-related Custom Fields
 Compound-related Custom Fields
 Custom Fields reported separately

Sequence Output
 Sequence Summary

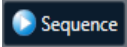

Select Destination for Printout

Printer
 File:

Path:
C:\Users\Public\Documents\ChemStation
\3\Sequence\

Exercise 2. Run the sequence

Now you are ready to acquire data with the sequence you just created.

Steps	Detailed Instructions	Comments
1 Confirm that your sequence includes four samples.	<ul style="list-style-type: none"> Verify that the Autosampler Tray diagram shows four samples. 	
2 Place the samples you prepared in Chapter 1 into the appropriate positions in the autosampler.		You prepared the samples in " Exercise 2. Prepare the samples for the analyses " on page 12.
3 Inject the samples.	<ul style="list-style-type: none"> Click Sequence in the Run Control Bar. 	This button is only available if you have selected Sequence mode on the main toolbar. 
4 (Optional) For the first blank analysis, monitor the total ion chromatogram and the UV chromatogram during data acquisition.	<ol style="list-style-type: none"> Activate the Online Plot window. Monitor the MS signal to ensure a stable baseline. 	As the sequence progresses, the Autosampler Tray diagram is color-coded as follows: Gray - samples that have been analyzed. White - samples not yet analyzed. Blue - the current sample.
5 When the sequence is done, view the Sequence Summary Report.	<ol style="list-style-type: none"> Retrieve the report from the printer. Examine the report to confirm that all the samples ran. 	
6 When the sequence is finished, view the results.	<ol style="list-style-type: none"> Display Data Analysis view. Load the first data file you just created. Examine the DAD and MS chromatograms. Repeat step b and step c for the other data files. 	<ul style="list-style-type: none"> If you need help, follow the general procedure in "Exercise 1. Display and manipulate chromatograms" on page 26 in Chapter 3. When you analyze your own samples, you can set up your method to automatically generate a data analysis report for each sample in the sequence.

This page intentionally left blank.

6

Quantitative Data Analysis

- Exercise 1. Create a method for quantification 58
 - Task 1. Create a new method 58
 - Task 2. Set up the signal for quantification 59
 - Task 3. Integrate the low-level standard 61
 - Task 4. Set general calibration parameters 63
 - Task 5. Set up the calibration curve 64
 - Task 6. Explore options to refine the calibration 68
- Exercise 2. Process a sample and print a report 69

This chapter shows you how to use the OpenLab ChemStation Data Analysis to perform quantification. The exercises in this chapter illustrate a simple calibration that uses data files that you received with your OpenLab ChemStation software.

Before you start


- Read the *Quick Start Guide*.
- Make sure that you have the caffeine data files on your OpenLab ChemStation. Check for the files under **C:\Users\Public\Documents\Chemstation\1\Data\MSDEMO**. The file names are **CafCal0x.D**, where **x** is a number from **1** to **5**.

Exercise 1. Create a method for quantification

In this exercise, you create a calibrated method that you can use to quantify caffeine in the demo data.

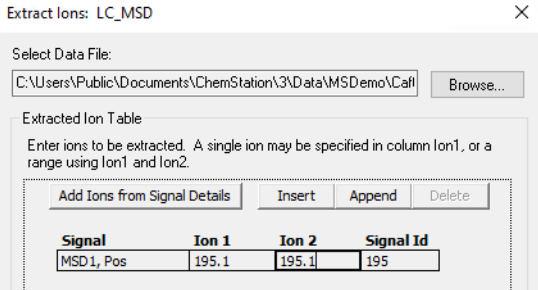

Task 1. Create a new method

In this task, you load a default method and save it to a new name. You later modify the new method to create a calibrated method.

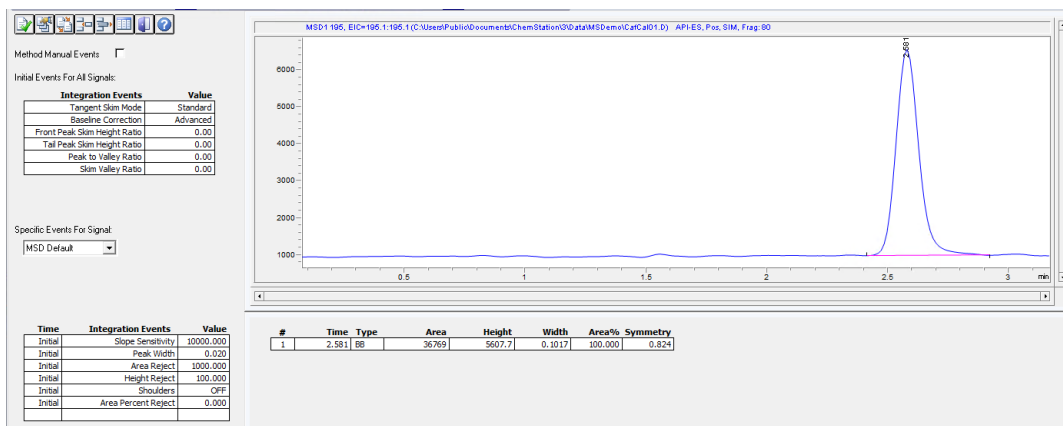
Steps	Detailed Instructions	Comments
1 Display Data Analysis view.	<ul style="list-style-type: none"> In the view selection area in the lower left of the OpenLab ChemStation window, click Data Analysis. 	
2 Load the method DEF_LC.M.	<ol style="list-style-type: none"> Click File > Load > Method. Navigate to the folder C:\Users\Public\Public Documents \Chemstation\1\Methods. Select the method file and click OK. 	
3 Save the method under the new name CAFFEINE CAL.M.	<ol style="list-style-type: none"> Select File > Save As > Method. Navigate to the folder C:\Users\Public\Public Documents \Chemstation\1\Methods. In the dialog box, for Name, type CAFFEINE CAL.M. Click OK. In the box for Comment for method history, type a comment. Click OK. 	

Task 2. Set up the signal for quantification

In this exercise, you add an extracted ion chromatogram (EIC) to the list of available signals for the method. Then you add this EIC to the Signal Details, so you can automatically load and integrate signals for the rest of the caffeine standards.

Steps	Detailed Instructions	Comments
1	<p>Open the data file CafCal01.D, located in the MSDEMO folder.</p> <p>a Select File > Load Signal.</p> <p>b Navigate to the folder: C:\Users\Public\Public Documents\Chemstation\1\Data\MSDEMO.</p> <p>c Select the data file CAFCAL01.D.</p> <p>d If necessary, clear the check box for Load using Signal Details.</p> <p>e In the Signals box, click the signal that begins with MSD1 TIC.</p> <p>f Click OK.</p>	
2	<p>Extract the major ion of caffeine.</p> <p>a Select File > Extract Ions.</p> <p>b For Ion 1, type 195.1.</p> <p>c For Ion 2, type 195.1.</p> <p>d Click OK.</p>	<ul style="list-style-type: none"> The 195 ion is the (M+H)⁺ ion.
		
3	<p>Display the Calibration Toolset.</p> <p>Click Calibration, which is near the center of the window.</p>	 Calibration

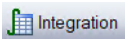



Steps	Detailed Instructions	Comments
4	<p>Set up the signal for quantification.</p> <p>a Do one of the following:</p> <ul style="list-style-type: none"> Click Edit to edit the current method signals. Select Calibration > Signal Details. <p>b From the list of Available Signals, select MSD1 195, EIC=195.1:195.1.</p> <p>c Click Add to Method.</p> <p>d Click OK.</p>	<p>The EIC signal is available only because you loaded the 195 EIC in step 2.</p>

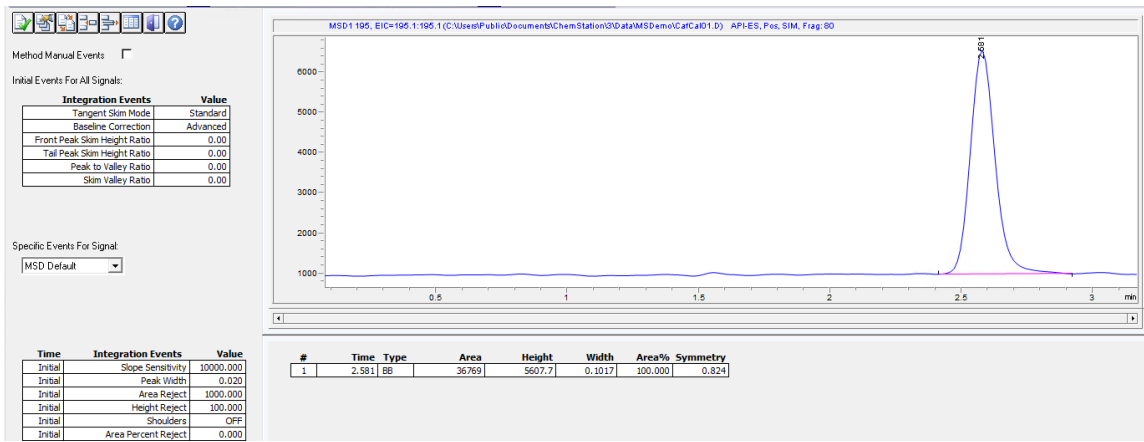



5	<p>(Optional) Save the method under the same name (CAFFEINE CAL.M).</p> <p>a Select File > Save > Method.</p> <p>b In the box for Comment for method history, type a comment.</p> <p>c Click OK.</p>	<p>For these exercises, you save the method frequently, but you could wait instead until you had established all the method settings.</p>
---	---	---

Task 3. Integrate the low-level standard

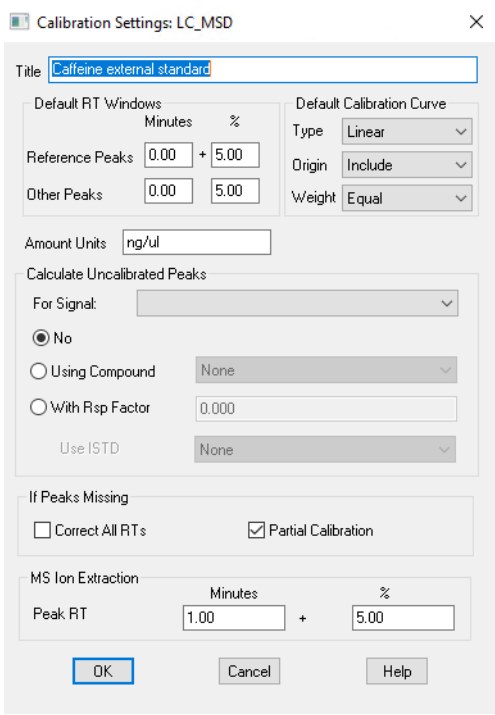
In this exercise, you establish integration parameters for your calibrated method. You use the low-level standard because it is usually the most difficult to integrate.

Steps	Detailed Instructions	Comments
1 Display the Integration Toolset.	<ul style="list-style-type: none"> Click Integration, which is near the center of the window. 	
2 Integrate the chromatogram.	<ul style="list-style-type: none"> a Click Auto Integrate, which is near the center of the window. b Check that you have five integrated peaks with these initial settings. 	<ul style="list-style-type: none"> Auto Integrate estimates initial integration parameters and then performs the integration.
3 Adjust the integration parameters to get one integrated peak.	<ul style="list-style-type: none"> a Click Edit/Set Integration Events Table. b In the integration events for all signals, for Baseline Correction, select Advanced. c Click Auto Integrate. d When you are prompted to save the events table, click Yes. e Verify that your results are the same or very similar to those shown below.  	<ul style="list-style-type: none"> For detailed information about integration events, see <i>OpenLab ChemStation Concepts and Workflows.pdf</i>.



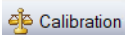

Steps	Detailed Instructions	Comments
4 Save the integration events with the method in memory.	<ul style="list-style-type: none">Click the button to exit and save the integration results.	
5 (Optional) Save the method under the same name (CAFFEINE CAL.M).	<ol style="list-style-type: none">Select File > Save > Method.In the box for Comment for method history, type a comment.Click OK.	

Task 4. Set general calibration parameters

Steps	Detailed Instructions	Comments
1	Establish calibration parameters.	<p>a Select Calibration > Calibration Settings.</p> <p>b In the Title box, type a title, for example <code>Caffeine external standard</code>.</p> <p>c Leave the rest of the items at the default settings, shown below.</p> <p>d Click OK.</p>
		
2	(Optional) Save the method under the same name (CAFFEINE CAL.M).	<p>a Select File > Save > Method.</p> <p>b In the box for Comment for method history, type a comment.</p> <p>c Click OK.</p>

Task 5. Set up the calibration curve

In this exercise, you integrate the rest of the standards and add all standards to the calibration curve.

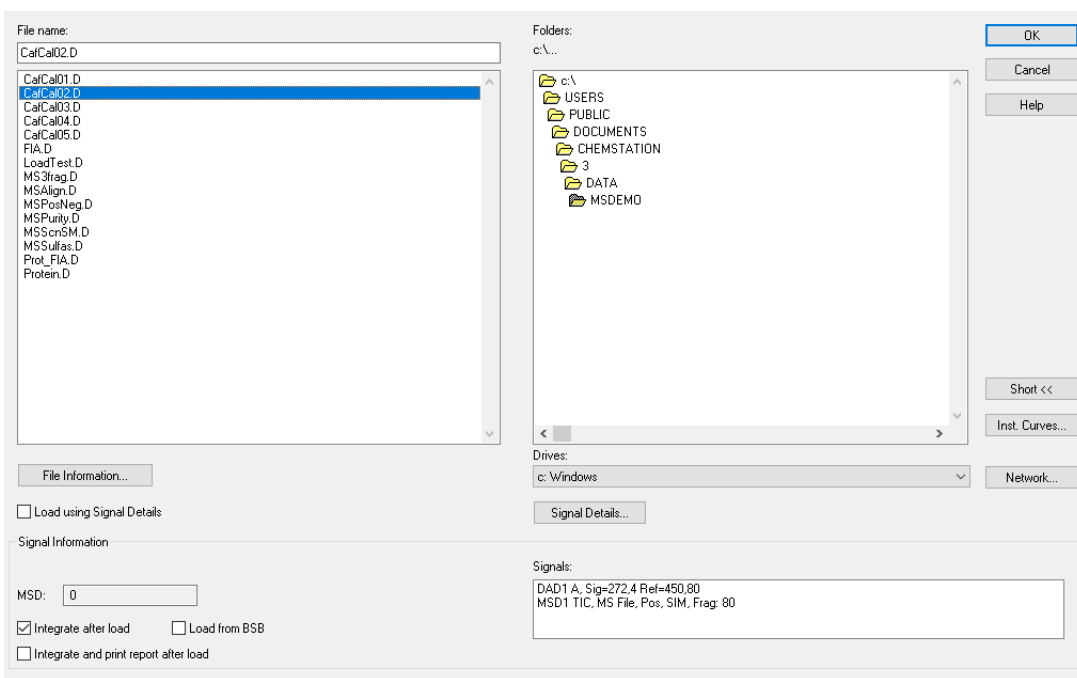
Steps	Detailed Instructions	Comments
1 Display the Calibration Toolset.	<ul style="list-style-type: none"> Click Calibration, which is near the center of the window. 	
2 Add the low-level standard to the calibration curve.	<p>a Do one of the following:</p> <ul style="list-style-type: none"> Click New Calibration Table. Select Calibration > New Calibration Table. <p>b Click Automatic Setup Level 1.</p> <p>c Click OK.</p> <p>d In the Calibration Table pane (shown below), under Compound, type caffeine and under Amt (amount), type 0.5.</p> 	<ul style="list-style-type: none"> Do not worry at this point if your calibration curve displays a message that says the curve is invalid.


Calibration Table

Enter Delete Insert... Print OK Help

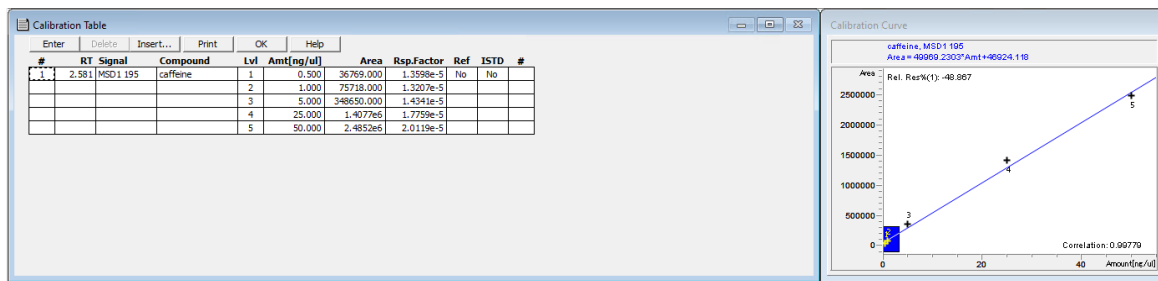
#	RT	Signal	Compound	Lvl	Amt[ng/ul]	Area	Rsp.Factor	Ref	ISTD	#
1	2.581	MSD1195	caffeine	1	0.500	36769.000	{ 1.3598e-5 }	No	No	

Steps	Detailed Instructions	Comments
3 Load and integrate the second standard.	<ol style="list-style-type: none"> Select File > Load Signal. Under File name, select CAFICAL02.D. Mark the check box for Load using Signal Details. Mark the check box for Integrate after load. Check that your dialog box looks like the one below. Click OK. 	<ul style="list-style-type: none"> These settings enable you in a single step to load the appropriate signal(s) and integrate them.



4 Add the second standard to the calibration curve.	<ol style="list-style-type: none"> Click Add new level. In the dialog box, for Default Amount, type 1 and click OK. Verify that the calibration table now has two entries, and the calibration curve contains two points. 	
---	---	---

Steps	Detailed Instructions	Comments
<p>5 Add the remaining three standards to the calibration table:</p> <ul style="list-style-type: none"> • CAFCAL03.D: 5 ng/μL • CAFCAL04.D: 25 ng/μL • CAFCAL05.D: 50 ng/μL 	<p>a Select File > Load Signal.</p> <p>b Under File name, select the next data file.</p> <p>c Verify that the chromatogram is properly integrated.</p> <p>d Click Add new level.</p> <p>e In the dialog box, for Default Amount, type the amount shown in step 5 and click OK.</p> <p>f Verify that the calibration table and the calibration curve contain the new entry.</p> <p>g Repeat step a through step f until you have added all the standards.</p> <p>h Confirm that your calibration table and calibration curve look like the ones below.</p>	<ul style="list-style-type: none"> • If multiple peaks are integrated in a chromatogram, retention time is used to find the correct peak for the calibration curve.



Steps	Detailed Instructions	Comments
6 Refine the calibration curve.	<p>a Select Calibration > Calibration Settings.</p> <p>b Under Default Calibration Curve, for Type, select Quadratic.</p> <p>c Click OK.</p> <p>d Verify that your calibration curve now looks like the one below.</p>	
	<p>The figure is a 'Calibration Curve' plot for 'caffeine, MSD1 195'. The y-axis is labeled 'Area' and ranges from 0 to 2,500,000. The x-axis is labeled 'Amount[ng/ul]' and ranges from 0 to 40. A blue quadratic curve is plotted through three data points labeled 3, 4, and 5. The equation for the curve is $Area = -263.12956 \cdot Amt^2 + 62452.865 \cdot Amt + 18663.094$. The correlation coefficient is 0.99991. The relative residual for the first point is -26.202.</p>	
7 (Optional) Save the method under the same name (CAFFEINE CAL.M).	<p>a Select File > Save > Method.</p> <p>b In the box for Comment for method history, type a comment.</p> <p>c Click OK.</p>	


Task 6. Explore options to refine the calibration

This exercise describes additional calibration table layouts that give you more calibration options. You do not need these options to process the caffeine demonstration data, but you may need them when you process your own samples.

Steps	Detailed Instructions	Comments
1 Explore options to change the way calibration curves are constructed.	<p>a Select Calibration Table Options > Peak Details.</p> <p>b Verify that you see these columns in the calibration table:</p> <ul style="list-style-type: none"> • Curve Type • Origin • Weight 	<ul style="list-style-type: none"> • Note that this calibration table layout lets you change: <ul style="list-style-type: none"> • Curve Type: The type of calibration curve (linear, quadratic, etc.) • Origin: How the origin (zero point) is treated. • Weight: The relative weights of the data points.
2 Explore options to add qualifier ions.	<p>a Select Calibration Table Options > Identification Details.</p> <p>b Verify that you see these columns in the calibration table:</p> <ul style="list-style-type: none"> • Resp % (response percent) • +- (window for the response percent) • Pk Usage (peak usage) 	<ul style="list-style-type: none"> • Note that this calibration table layout lets you define: <ul style="list-style-type: none"> • Pk Usage: How the calibration uses the peak, for example, as a main calibration ion or a qualifier ion • Resp %: The expected response of the qualifier ion, as a percentage of the main peak • +-: A window for the expected percentage.
3 Display the original options for the calibration table.	<p>a Select Calibration Table Options > Overview.</p> <p>b Verify that the calibration table looks the same as in step 5 on page 66.</p>	

Exercise 2. Process a sample and print a report

In this exercise, you specify a report and test your calibration method by processing one of the standards as if it were a sample. You print a report of the results.

Steps	Detailed Instructions	Comments	
1	<p>Specify a report with the following settings:</p> <ul style="list-style-type: none"> Report destination: Screen External standard (ESTD) calculation, based on area Report style: Detail 	<p>a Do one of the following:</p> <ul style="list-style-type: none"> Select Report > Specify Report. Click Specify Report. <p>b Enter parameters as described in step 1 and shown in the next two figures.</p> <p>c Click Quantitation settings.</p> <p>d Click OK.</p>	

Specify Report: LC_MSD

Reporting settings Quantitation settings

Use Intelligent Reporting Use Classic Reporting

Style

Report style:

Quantitative result sorted by:

Sample info on each page Add fraction table and ticks

Add chromatogram output Add summed peaks table

Add sample custom fields to sample info Add compound custom fields

Report layout for uncalibrated peaks

Separately With calibrated peaks Do not report

Chromatogram output

Portrait Landscape Multi-page

Size (% of page)

Time:

Response:

Pages

Destination

Printer Screen File

File setting

File prefix:

PDF CSV HTM DIF

Unique PDF file name TXT XLS EMF

Steps	Detailed Instructions	Comments						
<div style="border: 1px solid black; padding: 5px;"> <p>Specify Report: LC_MSD ×</p> <p>Reporting settings Quantitation settings</p> <div style="border: 1px solid gray; padding: 5px;"> <p>Calculation mode</p> <p>Calculate: ESTD Based on: Area</p> </div> <div style="border: 1px solid gray; padding: 5px; margin-top: 5px;"> <p>ISTD Correction</p> <p><input checked="" type="checkbox"/> Use multiplier and dilution factor with ISTDs</p> </div> <div style="border: 1px solid gray; padding: 5px; margin-top: 5px;"> <p>Calculation factors</p> <p>Use sample data: From data file</p> <p>Sample amount: 0</p> <p>Multiplier: 1</p> <p>Dilution: 1</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th>#</th> <th>Compound</th> <th>ISTD Amount</th> </tr> </thead> <tbody> <tr style="background-color: #cccccc;"> <td colspan="3" style="height: 40px;"> </td> </tr> </tbody> </table> </div> <div style="text-align: right; margin-top: 10px;"> OK Cancel Help </div> </div>			#	Compound	ISTD Amount			
#	Compound	ISTD Amount						

- 2 Save the method under the same name (CAFFEINE CAL.M).
 - a Select **File > Save > Method**.
 - b In the box for **Comment for method history**, type a comment.
 - c Click **OK**.

- 3 Load the standard of medium concentration.
 - a Select **File > Load Signal**.
 - b Under **File name**, select **CAFAL03.D**.

Steps	Detailed Instructions	Comments
4 Process the medium-level standard and print the report.	<p>a Do one of the following:</p> <ul style="list-style-type: none"> • Select Report > Print Report. • Click preview results. <p>b Verify that page 1 of the report contains header information, an integrated chromatogram, and an external standard report.</p> <p>c Check that the caffeine amount is about 5 ng/μL.</p> <p>d At the bottom of the report window, click Next.</p> <p>e Verify that page 2 of the report shows the calibration curve with the measured point identified with dotted lines.</p> <p>f (Optional) At the bottom of the report window, click Print so you get a hard copy.</p> <p>g At the bottom of the report window, click Close.</p>	<ul style="list-style-type: none"> • Another way to generate a hard copy is to click Print Report.



In This Book

When you do the exercises in this book, you learn how to:

- Prepare your MSD system for an analysis
- Set up methods for scan and selected ion monitoring analyses
- Acquire data
- Set up sequences for automated sample analyses
- Perform qualitative and quantitative analyses.

www.agilent.com

© Agilent Technologies, Inc. 2022

June 2022



D0015483

