

Agilent MassHunter BioConfirm Software

Familiarization Guide

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

Basic Tasks 3

- Task 1. Open the BioConfirm program 3
- Task 2. Zoom in and out of the chromatogram 6
- Task 3. Change window layouts 8
- Task 4. Creating a Protein Sequence File 10

Intact Protein Workflow 12

- Exercise 1. Interactive Intact Protein Workflow 13
- Exercise 2. Automated Intact Protein Workflow 17

Protein Digest Workflow 19

- Exercise 3. Interactive Protein Digest Sequence Matching 20
- Exercise 4. Automated Protein Digest Workflow 25

Released Glycans Workflow 26

- Exercise 5. Interactive Released Glycans 27
- Exercise 6. Automated Released Glycans Workflow 29

Review Results 33

- Exercise 7. Reprocessing Samples 33
- Exercise 8. Using Result Review mode 36
- Exercise 9. Using Report Builder 40

Deconvolution 45

- Exercise 10. Interactive Protein Molecular Weight Determination 45
- Exercise 11. Using the Mirror Plot window 49
- Exercise 12. Viewing Biomolecule Information 53

Where to find more information

- *Agilent MassHunter BioConfirm Software Quick Start Guide*
- Agilent MassHunter BioConfirm eFamiliarization
- Agilent MassHunter BioConfirm Training Videos
- Online Help provides in-depth information and can be displayed in the following ways:
 - Click **Contents** or **Search** from the BioConfirm software Help menu.
 - Press the **F1** key to get more information about a window or dialog box.

How to use this guide

Try to do these familiarization exercises initially using the steps listed in the first column. Then if you need more information, follow the detailed instructions in the second column.

Before you start

Copy the data files used for these tasks onto your hard disk as follows:

- 1 Copy all of the data files from the **Data** folder on the BioConfirm setup media to your computer hard drive. We recommend copying the data files to the **D:\MassHunter\Data** folder.
- 2 Make sure you have both read and write permissions for the folder you just created on your computer. This is required if you want to save results.
 - a In Windows Explorer right-click the folder where you copied the data files and click **Properties** from the shortcut menu.
 - b *Clear* the **Read-only Attributes** check box if it is marked.
 - c In the Confirm Attribute Changes dialog, click **Apply changes to this folder, subfolders, and files**, and then click **OK**.
- 3 Copy all of the sequences from the **ProteinSequences** folder on the BioConfirm setup media to your computer hard drive. We recommend copying the data files to the **D:\MassHunter\ProteinSequences** folder.

Basic Tasks

Task 1. Open the BioConfirm program

Basic Tasks

Task 1. Open the BioConfirm program

In this task you open multiple data files using the current method.

Task 1. Open the BioConfirm program with multiple data files

Steps	Detailed Instructions	Comments
<p>1 Open the BioConfirm program.</p> <ul style="list-style-type: none">Open these data files NIST mAb 1.d Nist mAb 2.d NIST mAb Digest.d NIST mAb Digest2.d ReleasedGlycans1.d ReleasedGlycans2.d in the folder D:\MassHunter\Data, or in the folder where you copied them.Make sure that the Use current method button is clicked.Make sure that the Load result data check box is cleared.	<p>a Double-click the Agilent MassHunter BioConfirm 10.0 icon. The system displays the Open Sample dialog box.</p> <p>b Go to the folder D:\MassHunter\Data or to the folder where the example files are located.</p>	<ul style="list-style-type: none">You can get help for any window, dialog box, or tab by pressing the F1 key when that window is active.

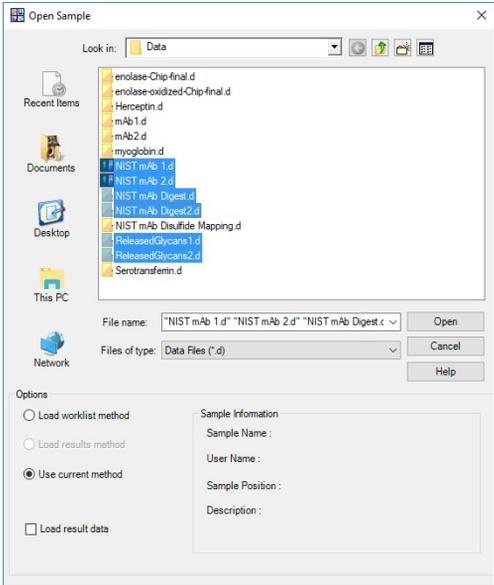


Figure 1. Open data files when opening software

Basic Tasks

Task 1. Open the BioConfirm program

Task 1. Open the BioConfirm program with multiple data files (continued)

Steps	Detailed Instructions	Comments
	<p>c Click the NIST mAb 1.d file.</p> <p>d Press and hold the Shift key while you click NIST mAb Digest.d.</p> <p>e Press and hold the Ctrl key while you click ReleasedGlycans1.d and ReleasedGlycans2.d.</p> <p>f Clear the Load result data check box.</p> <p>g Click Open. All the data files are displayed in the Sample Table window. The selected sample in the Sample Table is also shown in the Sample Chromatogram Results window.</p> <p>h Click the List Mode button in the Sample Chromatogram Results toolbar.</p> <p>i Click the NIST mAb 1.d data file.</p>	<ul style="list-style-type: none">• If you press the Shift key, you can pick a group of files that are directly next to each other.• If you press the Ctrl key, you can pick files which are not directly next to each other in the list.• What you see in the main window at this point depends on the method, layout, display and plot settings used before you opened these files.• When you click the List Mode button, the background of the button changes to orange.
2 Return the main window to the default Intact Protein layout.	<ul style="list-style-type: none">• Click Intact Protein Layout in the main toolbar.	<ul style="list-style-type: none">• You click the  button in the graphics window to change the display options.• You can switch between layouts for the different workflows when you click the buttons in the main toolbar.• You can change the layout if you click Configuration > Window Layouts > Load Layout.

Basic Tasks

Task 1. Open the BioConfirm program

Task 1. Open the BioConfirm program with multiple data files (continued)

Steps

Detailed Instructions

Comments

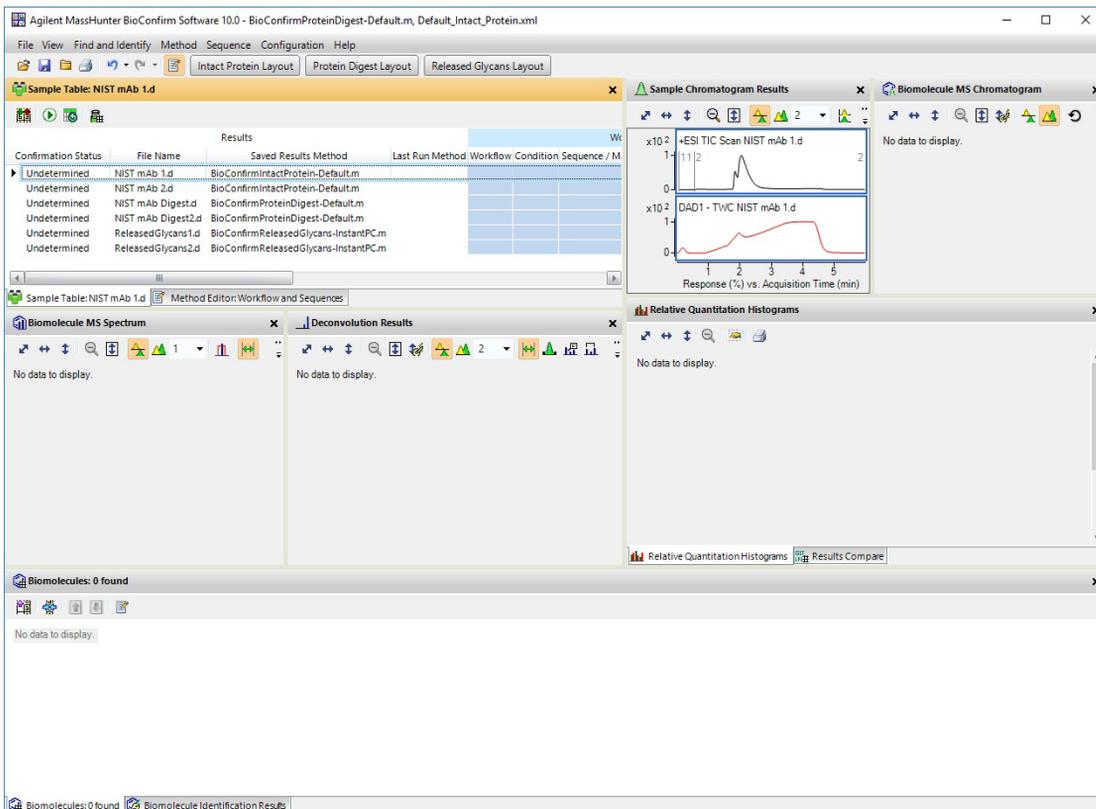


Figure 2. BioConfirm main window

Basic Tasks

Task 2. Zoom in and out of the chromatogram

Task 2. Zoom in and out of the chromatogram

In this task, you become familiar with the zoom in and zoom out features of the BioConfirm program.

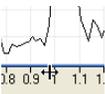
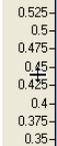
Task 2. Zoom in and out of the chromatogram

Steps	Detailed Instructions	Comments
1 Practice zooming in and out on the chromatogram in the Sample Chromatogram Results window. <ul style="list-style-type: none">• Zoom in twice on the peak.• Zoom in one more time autoscaling the y-axis.• Zoom out once to the previous zoom position.• Completely zoom out to the original chromatogram.	<p>a Click the right mouse button and drag over an area on the last peak. Make sure that the Autoscale Y-axis during Zoom button, , is not selected for this step.</p> <p>b Repeat step a.</p> <p>c Click the Autoscale Y-axis during Zoom button, , in the toolbar.</p> <p>d Click the right mouse button again and drag over an area of the peak for the third time. The BioConfirm program automatically scales the y-axis to the largest point in the range. </p> <p>e Click the Unzoom button, , to undo the last zoom operation. You can undo the last fifteen zoom operations.</p> <p>f Click the Autoscale X-axis and Y-axis button, , to zoom out completely.</p>	<ul style="list-style-type: none">• You can also use these zoom features in the Biomolecule MS Spectrum window, the Biomolecule Fragment Spectrum window, the Deconvolution Results window, the Deconvolution Mirror Plot window, and the Biomolecule MS Chromatogram window.• In addition to those windows, you can also zoom on the x-axis and y-axis and use the toolbar buttons in the Relative Quantitation Histograms window. You cannot drag over an area in the Relative Quantitation Histograms window.• A selected button has an orange background color.

Basic Tasks

Task 2. Zoom in and out of the chromatogram

Task 2. Zoom in and out of the chromatogram (continued)

Steps	Detailed Instructions	Comments
<p>2 Practice zooming in and out on each axis separately.</p> <ul style="list-style-type: none">Zoom in only along the x-axis. <p>Hint: Right-click the x-axis values and move cursor from left to right.</p> <ul style="list-style-type: none">Partially zoom out the x-axis. <p>Hint: Move cursor in opposite direction.</p> <ul style="list-style-type: none">Completely zoom out of the x-axis.Repeat the previous steps for the y-axis.	<p>a To zoom in on the x-axis, move the cursor to the x-axis values until a horizontal double arrow appears.</p> <p>b Click the right mouse button and drag the new cursor from left to right across the x-axis values.</p> <p>c To zoom out on the x-axis, click the right mouse button and drag from right to left on the x-axis values.</p> <p>d Click the Autoscale X-axis button, , to completely zoom out on the x-axis.</p> <p>a To zoom in on the y-axis, move the cursor to the y-axis values until a vertical double arrow appears.</p> <p>b Click the right mouse button and drag the new cursor from bottom to top across the y-axis values.</p> <p>c To zoom out on the y-axis, click the right mouse button and drag from the top towards the bottom of the y-axis values.</p> <p>d Click the Autoscale Y-axis button, , to completely zoom out on the y-axis.</p>	<p> Horizontal Double Arrow</p> <p> New cursor appears when you right-click the x-axis value</p> <p> Vertical Double Arrow</p> <p> New cursor appears when you right-click the y-axis values.</p>

Basic Tasks

Task 3. Change window layouts

Task 3. Change window layouts

In this task, you move windows within the main view and create various window layouts. Default layouts are available for each workflow.

Task 4. Change window layout

Steps	Detailed Instructions	Comments
<p>1 Change the window layout:</p> <ul style="list-style-type: none">• Change the window size.• Save a window layout.• Unlock the layout.• Change the Chromatogram Results window to be floating.• Move the Chromatogram Results window.• Display the tools for repositioning the windows.	<ul style="list-style-type: none">• To change the size of a window, drag the boundary between the windows.• To load the default layout for a workflow, click one of the buttons in the main toolbar: Intact Protein Layout, Protein Digest Layout, and Released Glycans Layout.• To load a layout, click Configuration > Windows Layouts > Load Layout.• To save a window layout, click Configuration > Window Layouts > Save Layout.• To lock or unlock a layout, click Configuration > Window Layouts > Lock Layout.• To make a window float, right-click the title bar of the window, and click Floating from the shortcut menu.• To move a window, click the title bar of the window and drag the window to the desired location.• To display the repositioning tools, drag the window over one of the other windows. When one window is overlapped with another, the program displays several layout tools, as shown in Figure 3.	<ul style="list-style-type: none">• If the layout is locked, the system displays a check mark next to the Lock Layout menu.• You can only use the repositioning tools when the layout is unlocked.• You can also make a window float by double-clicking the title bar of the window.• The following layouts are shipped with the software: Default_IntactProtein.xml Default_Protein_Digest.xml Default_Released_Glycans.xml

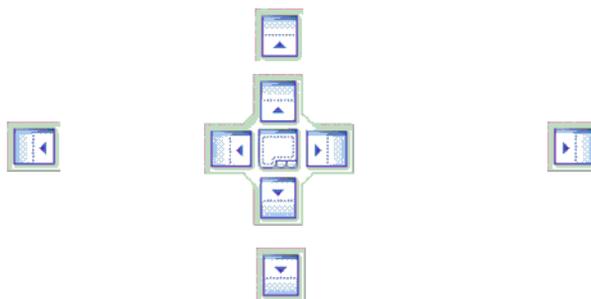


Figure 3. Window repositioning tools

Basic Tasks

Task 3. Change window layouts

Task 4. Change window layout (continued)

Steps	Detailed Instructions	Comments
<p>2 Reposition the Sample Chromatogram Results window.</p> <ul style="list-style-type: none">• Move the window so that it is at the top, to the left, to the right and then at the bottom of the other windows.• Move two windows together so that they are on top of one another and available only through the tabs at the bottom.• Restore the default layout.	<ul style="list-style-type: none">• If you drag the cursor over one of the smaller icons, the window you are dragging will be placed above, to the right, below, or to the left of all of the other windows.• Drag the cursor over the larger icon. The window can also be placed above, to the right, below, or to the left of the other window by dragging the cursor over the edges of the larger icon.• To tab two windows together, drag the cursor over the center of the larger icon. You will see a shadow version of the two windows tabbed together. Stop dragging the mouse. The two windows will be tabbed together.• Click Intact Protein Layout in the main toolbar.	<ul style="list-style-type: none">• The cursor must be over one of the arrows in a box in order for repositioning to occur.• Clicking the Configuration > Load Default Layout command restores the default layout. The default layout is different than the default layout for each workflow.

Basic Tasks

Task 4. Creating a Protein Sequence File

Task 4. Creating a Protein Sequence File

This task guides you through the creation of a myoglobin sequence file.

Steps	Detailed Instructions	Comments
1 Start the Agilent MassHunter Sequence Manager.	<ul style="list-style-type: none">Click Sequence > Sequence Manager.	
2 Create a new sequence.	<ul style="list-style-type: none">a Type <code>Myoglobin</code> for the name of the Sequence.b Click the + button. The Sequence Editor pane opens automatically with a new sequence displayed for editing.	<ul style="list-style-type: none">Protein is automatically selected for the sequence type.
3 Enter the amino acid sequence shown below into the Sequence Manager.	<ul style="list-style-type: none">Type in individual amino acids one at a time between the N-term and C-term symbols. <p>GLSDGEWQQVLNVWVGKVEADIAGHGQEVLRIRLFTGHPETLEKFDKFKHLKTEAEM KASEDLKKHGTVVLTAALGGILKKKGHHEAELKPLAQSHATKHKIPIKYLEFISDAIIH VLHSHKHPGDFGADAQGAMTKALELFRNDIAAKYKELGFQG</p>	<ul style="list-style-type: none">Use the single-character (letter) amino acids abbreviations.Tip: If you are reading this document as a PDF file on your computer, you can copy and paste the sequence into the Sequence Manager window.

Note: The myoglobin sequence does not have any links or modifications, but some sequences do. In that case, add links and modifications as described in the *Quick Start Guide* or *online Help*.

Basic Tasks

Task 4. Creating a Protein Sequence File

Steps	Detailed Instructions	Comments
4 Save the sequence as the name <i>iii_myoglob.psq</i> , where <i>iii</i> represents your initials.	<p>a Click Sequence > Export Sequences.</p> <p>b Type <i>iii_myoglob</i> in the File name box.</p> <p>c Click Save.</p>	<ul style="list-style-type: none">The sequence is saved as a .psq file that can be imported for use in other methods as described in Exercise 4 or referenced from worklists as described in Exercise 5.

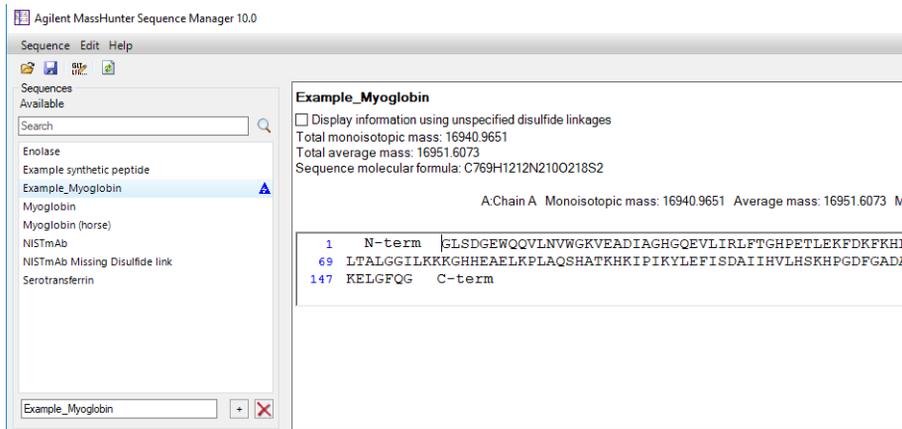


Figure 4. Creating a sequence file of myoglobin in the Sequence Manager program

Intact Protein Workflow

Task 4. Creating a Protein Sequence File

Intact Protein Workflow

Step 1 - Open the data file of interest.

Step 2 - Open a BioConfirm Method or create a new one.

Step 3 - Edit sequences if necessary in the Sequence Manager program:

- Add or edit the sequence text.
- Apply or edit modifications
- Apply or edit links

Step 4 - Select **Intact Protein** for the **Workflow** on the **Workflow and Sequences** tab. Select the **Condition**.

Step 5 - Select the **Sequence/Masses** to match on the Workflow and Sequences tab.

If the sequence you want to match is not in the method or Sample Table, then:
Import or create a sequence.

Step 6 - Select the **Mods and Profiles** on the Workflow and Sequences tab.

Step 7 - Run the Method Workflow.

Step 8 - Review the results which are shown in these windows:

Sample table

Biomolecules table

Biomolecule Identification Results

Deconvolution Results

Biomolecule MS Chromatogram

Biomolecule MS Spectrum

Biomolecule Fragment Spectrum

Results Compare

Relative Quantitation Histograms

Step 9 - Print report.

Exercise 1. Interactive Intact Protein Workflow

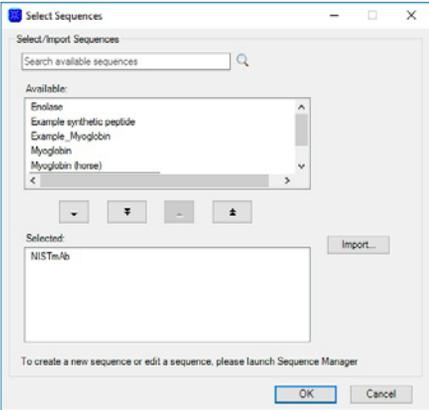
This exercise shows you how to set method parameters, match an intact protein sequence, and view the results. This exercise uses the **NISTmAb.seq** sequence file and the **NIST mAb1.d** data file copied before you started. See **“Before you start”** on page 2.

If you select the Intact Protein workflow, the Find by Protein Deconvolution algorithm runs and uses protein Matching Rules (Intact Protein, and Predicted Modifications). You can select whether or not Protein Truncation is done.

Steps	Detailed Instructions	Comments
1 Open the method to use as a starting point for the new method.	<ol style="list-style-type: none"> a Click Method > Open. b Select the BioConfirmIntactProtein-Default.m folder. c Click Open. 	
2 If the NIST mAb1.d data file is not already open, open it.	<ol style="list-style-type: none"> a Click File > Open Data File. b Locate the NIST mAb1.d folder. c Click Open. 	<ul style="list-style-type: none"> • The TIC is automatically displayed in the Sample Chromatogram Results window.
3 Display the Deconvolute (Protein) section in the Method Editor window.	<ol style="list-style-type: none"> a Click View > Method Editor if the Method Editor is not visible. b Select Intact Protein > Deconvolute (Protein) in the Method Editor window. 	
4 Run the Find by Protein Deconvolution algorithm.	<ol style="list-style-type: none"> a Review the settings and modify them if necessary. b Click  on the Method Editor toolbar to start the Find by Protein Deconvolution algorithm. c If the Find Proteins dialog box opens, select NIST mAb 1.d and click Find. d Review the results in the Biomolecules window. 	<ul style="list-style-type: none"> • In this case you are using the default method parameters. For some data files, you will need to use different parameters as described in the <i>Quick Start Guide</i> or <i>online Help</i>. • If you have more than one data file open, the Find Proteins dialog box opens.
5 Display the Workflow and Sequences section in the Method Editor window.	<ul style="list-style-type: none"> • Click Method Automation > Workflow and Sequences in the Method Editor window. 	

Intact Protein Workflow

Exercise 1. Interactive Intact Protein Workflow

Steps	Detailed Instructions	Comments
6 Import the NISTmAb sequence.	<ol style="list-style-type: none">Select Intact Protein for the Workflow.Select non-reduced for the Condition.Click the  button next to the Sequences parameter. The Select Sequences dialog box opens.Double-click NISTmAb. If NISTmAb is not available, click Import.Select NISTmAb.psq and click Open.Verify that the NISTmAb sequence is in the Selected list.Click OK.	<ul style="list-style-type: none">You will use the sequence as is. You can add or modify modifications and links to sequences as described in <i>online Help</i> and the <i>Quick Start Guide</i>.
		
7 Select the mAb modification.	<ol style="list-style-type: none">Click the  button next to the Modifications parameter. The Select Modifications and Profiles dialog box opens.Double-click mAb in the Available list in the Modifications and profiles section.Click OK.	<ul style="list-style-type: none">The mAb sequence has modifications. You can learn how to add modifications in the <i>online Help</i> and the <i>Quick Start Guide</i>.
8 Start the match search.	<ol style="list-style-type: none">Click Intact Protein > Match Tolerances.Click  on the Method Editor toolbar.Select NIST mAb 1.d and click Match.	<p>Alternate methods:</p> <ul style="list-style-type: none">Click Find and Identify > Match Sequences.
9 Review the results.	<ul style="list-style-type: none">Select the Biomolecule 1 row in the Biomolecules table.	<ul style="list-style-type: none">In the BioConfirmIntactProtein-default layout, the Biomolecule Identification Results window is tabbed with the Biomolecules window.

Intact Protein Workflow

Exercise 1. Interactive Intact Protein Workflow

Steps

Detailed Instructions

Comments

The screenshot displays the Agilent MassHunter BioConfirm Software 10.0 interface. The main window is titled "BioConfirmIntactProtein-Default.m, Default_Intact_Protein.xml". The "Sample Table" shows a list of results for "NIST mAb 1.d", including "Confirmed", "Undetermined", and "Released Glycans" entries. The "Sample Chromatogram Results" panel shows two chromatograms: "+ESI TIC Scan NIST mAb 1.d" and "DAD1 - TWC NIST mAb 1.d". The "Biomolecule MS Chromatogram" shows a peak at 4.611 minutes. The "Biomolecule MS Spectrum" shows a mass spectrum with peaks at 1538.9916, 1483.301, 2906.9171, and 4841.5318. The "Deconvolution Results" panel shows two deconvoluted mass spectra. The "Results Compare" panel shows a table of results for "NISTmAb" with columns for "Sequence/Mass", "Confirmation Status", "File", "RT (Range)", "Label", "Peak Mass (Da)", and "Pred M". The "Biomolecules: 32 found" table is shown below, with columns for "Label", "Mass", "RT", "Height", "Area", "Score", "Min Z", "Max Z", "ID Techniques Applied", "File", "ID Source", "Mining Algorithm", "Use for %Quant", "Area (MS)", and "%Quant".

Label	Mass	RT	Height	Area	Score	Min Z	Max Z	ID Techniques Applied	File	ID Source	Mining Algorithm	Use for %Quant	Area (MS)	%Quant
Biomolecule 1: A(1482013347	2.025	2488	9166272	92.13	30	101	Sequence Match	NIST mAb 1.d	Sequence Match	Maximum Entropy Deconv	<input checked="" type="checkbox"/>	1765983	11.26
Biomolecule 2: A(1483633865	1.859	2335	1336501	85.49	30	90	Sequence Match	NIST mAb 1.d	Sequence Match	Maximum Entropy Deconv	<input checked="" type="checkbox"/>	1737180	11.07
Biomolecule 3: A(1480393039	1.842	1567	904188	97.74	30	81	Sequence Match	NIST mAb 1.d	Sequence Match	Maximum Entropy Deconv	<input checked="" type="checkbox"/>	1039079	6.62
Biomolecule 4: A(148524599	1.842	1428	1034340	48.52	30	69	Sequence Match	NIST mAb 1.d	Sequence Match	Maximum Entropy Deconv	<input checked="" type="checkbox"/>	1266461	8.07
Biomolecule 5: A(14832409	2.107	1138	211834	0	118	185	Sequence Match	NIST mAb 1.d	Sequence Match	Maximum Entropy Deconv	<input checked="" type="checkbox"/>	1290098	8.22
Biomolecule 6	1481605959	2.058	1078	1792174	0	30	106		NIST mAb 1.d		Maximum Entropy Deconv	<input type="checkbox"/>	868267	
Biomolecule 7	1484858489	1.859	1024	738156	0	30	74		NIST mAb 1.d		Maximum Entropy Deconv	<input type="checkbox"/>	707014	

10 Save the method for use in Exercise 2.

- Click **Method > Save As**.
- Type the **File name** *iii_NIST_mAb.m*, where *iii* represents your initials.
- Click **Save**.

11 Investigate the Relative Quantitation feature. You normally use this feature to quantitate proteoforms (PTMs on the protein).

- In the Biomolecules window, hide all empty columns. Click  in the Biomolecules toolbar.
 - Review the values for **Use for %Quant** column for the biomolecules.
 - Review the values for the **%Quant (Height)** and **%Quant (Area)**.
- You can right-click the window and click **Add/Remove** columns to change the columns that are available.

Exercise 2. Automated Intact Protein Workflow

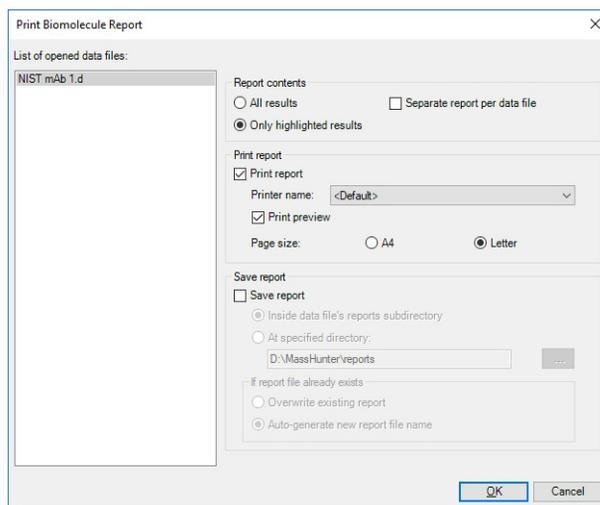
This exercise guides you through the setup of a worklist to automatically confirm the presence of NISTmAb in a previously acquired sample. This exercise uses the **NIST mAb1.d** data file copied in Exercise 1.

Steps	Detailed Instructions	Comments
1 If not already open, open the method <i>iii_NIST_mAb_Intact.m</i> .	<ul style="list-style-type: none"> a Click Method > Open. b Select the <i>iii_NIST_mAb_Intact.m</i> folder. c Click Open. 	This method was created in “ Exercise 1. Interactive Intact Protein Workflow ” on page 13.
2 Open the automation section in the Method Editor window.	<ul style="list-style-type: none"> • Click Method Automation > Workflow and Sequences in the Method Editor window. 	
3 Use the Intact Protein Workflow.	<ul style="list-style-type: none"> • Confirm that Intact Protein is selected for the Workflow. 	<ul style="list-style-type: none"> • In this case you are using the default method parameters. For some data files, you will need to use different parameters as described in the <i>Quick Start Guide</i> or <i>online Help</i>.
4 Import the NISTmAb sequence.	<ul style="list-style-type: none"> a Select non-reduced for the Condition. b Click the  button next to the Sequences parameter. The Select Sequences dialog box opens. c If NISTmAb is not available, click Import. d Select <i>NISTmAb.psq</i> and click Open. e Verify that the NISTmAb sequence is in the Selected list. f Click OK. 	<ul style="list-style-type: none"> • The NISTmAb.psq sequence file is available on the BioConfirm setup media. • You can learn about modifications and links in the <i>online Help</i> and in the <i>Quick Start Guide</i>.
5 Save the method.	<ul style="list-style-type: none"> • Click Method > Save. 	
6 Run the method workflow or run method automation.	<ul style="list-style-type: none"> • Click Method > Run Method Workflow. • Click Method > Run Method Automation (Workflow + Reports). • Click  on the Method Editor toolbar. • Reprocess the sample. See “Exercise 7. Reprocessing Samples” on page 33. 	<ul style="list-style-type: none"> • Method Automation first runs the method workflow, and then extracts additional chromatograms and generates a biomolecule report and exports results.

Intact Protein Workflow

Exercise 2. Automated Intact Protein Workflow

Steps	Detailed Instructions	Comments
7 (optional) Review the printed Biomolecule reports.	<ul style="list-style-type: none">If you clicked Run Method Automation, then a report is generated automatically.You can click File > Print > Biomolecule Report to generate a report for the current sample.	<ul style="list-style-type: none">You set report options in the Method Editor window in the Method Automation > Reports section.If you are creating a report interactively, you specify some of these parameters in the Print Biomolecule Report dialog box.



Protein Digest Workflow

Exercise 2. Automated Intact Protein Workflow

Protein Digest Workflow

The steps outlined below show the workflow for Protein Digest with Agilent MassHunter BioConfirm Software.

Step 1 - Open the data file of interest.

Step 2 - Open a BioConfirm Method or create a new one.

Step 3 - Edit sequences if necessary in the Sequence Manager program:

- Add or edit the sequence text.
- Apply or edit modifications
- Apply or edit links

Step 4 - Select the **Workflow** on the **Workflow and Sequences** tab. Select the **Condition**.

Step 5 - Select the **Sequence/Masses** to match on the Workflow and Sequences tab.

If the sequence you want to match is not in the method or Sample Table, then: Import or create a sequence.

Step 6 - Select the **Mods and Profiles** on the Workflow and Sequences tab.

Step 7 - Mark the **Enzymes** on the Workflow and Sequences tab.

Step 8 - Run the Method Workflow.

Step 8 - Review the results which are shown in these windows:

Biomolecules table
Biomolecule Identification Results
Sequence Coverage Map
Biomolecule MS Spectrum
Biomolecule Fragment Spectrum
Peptide Relative Quantitation Results
Results Compare
Relative Quantitation Histograms

Step 9 - Print report.

Exercise 3. Interactive Protein Digest Sequence Matching

This exercise shows you how to confirm protein digests interactively.

If you select the Protein Digest workflow, the Find Peptides algorithm runs and uses the enzyme selected in the Workflow and Peptides section and then runs the Protein Digest matching rules. See **“Before you start”** on page 2.

Steps	Detailed Instructions	Comments
1 Open the method to use as a starting point for the new method.	<ol style="list-style-type: none"> a Click Method > Open. b Select the BioConfirmProteinDigest-Default.m folder. c Click Open. 	<ul style="list-style-type: none"> • The parameters in the BioConfirmProteinDigest-Default.m method are a good starting point for Protein Digests.
2 Load the Protein Digest default layout.	<ul style="list-style-type: none"> • Click Protein Digest Layout on the main toolbar. 	<ul style="list-style-type: none"> • The TIC is automatically displayed in the Sample Chromatogram Results window.
3 Select NIST mAb Digest.d. If necessary, open the example sample file.	<ul style="list-style-type: none"> • If available, select NIST mAb Digest.d. otherwise a Click File > Open Data File. b Locate the NIST mAb Digest.d sample. c Clear the Load result data check box. d Click Open. 	<ul style="list-style-type: none"> • The TIC is automatically displayed in the Sample Chromatogram Results window.
4 Review the parameters in the Find Peptides section in the Method Editor window.	<ol style="list-style-type: none"> a Select Protein Digest > Find Peptides in the Method Editor window. b Review the settings on the various tabs of the Find Peptides section. c Click the MS-Only Extraction tab. d Review the parameters. For the example file, you can restrict the mass range to 300 - 1700. e In the MS-Only Extraction tab, enter 500 for the Use peaks with height >= counts. 	<ul style="list-style-type: none"> • You can change the default parameters as described in the next steps. You can also use the method without any changes. • For some data files, you will need to use different parameters as described in the <i>online Help</i>. • A very low peak height filter can result in greater sequence coverage but requires much more time to process.
5 Find biomolecules.	<ol style="list-style-type: none"> a Click  on the Method Editor toolbar to start the biomolecule search. b If the Find Peptides dialog box opens, select NIST mAb Digest.d and click Find. c When processing is complete, review the results in the Biomolecules window. 	<ul style="list-style-type: none"> • You can instead click Find and Identify > Find Peptides.

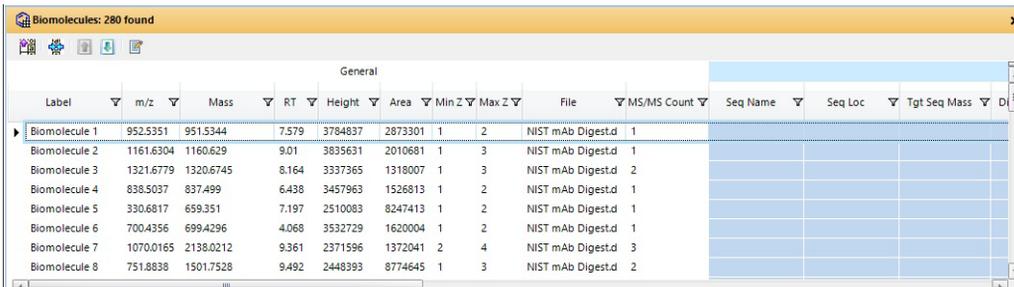
Protein Digest Workflow

Exercise 3. Interactive Protein Digest Sequence Matching

Steps

Detailed Instructions

Comments



The screenshot shows a software window titled "Biomolecules: 280 found". It contains a table with columns: Label, m/z, Mass, RT, Height, Area, Min Z, Max Z, File, MS/MS Count, Seq Name, Seq Loc, Tgt Seq Mass, and Di. The table lists 8 biomolecules with their respective values.

Label	m/z	Mass	RT	Height	Area	Min Z	Max Z	File	MS/MS Count	Seq Name	Seq Loc	Tgt Seq Mass	Di
Biomolecule 1	952.5351	951.5344	7.579	3784837	2873301	1	2	NIST mAb Digest.d	1				
Biomolecule 2	1161.6304	1160.629	9.01	3835631	2010681	1	3	NIST mAb Digest.d	1				
Biomolecule 3	1321.6779	1320.6745	8.164	3337365	1318007	1	3	NIST mAb Digest.d	2				
Biomolecule 4	838.5037	837.499	6.438	3457963	1526813	1	2	NIST mAb Digest.d	1				
Biomolecule 5	330.6817	659.351	7.197	2510083	8247413	1	2	NIST mAb Digest.d	1				
Biomolecule 6	700.4356	699.4296	4.068	3532729	1620004	1	2	NIST mAb Digest.d	1				
Biomolecule 7	1070.0165	2138.0212	9.361	2371596	1372041	2	4	NIST mAb Digest.d	3				
Biomolecule 8	751.8838	1501.7528	9.492	2448393	8774645	1	3	NIST mAb Digest.d	2				

6 Import the sequence.

- Click **Method Automation > Workflow and Sequences** in the Method Editor window.
- Select **Protein Digest** as the **Workflow**.
- Select **reduced** as the **Condition**.
- Click the next to the **Sequences/Masses** parameter.
- Double-click **NISTmAb**.
- Click **OK** in the **Select Sequences** dialog box.
- Click the next to the **Mods and Profiles** parameter.
- Double-click **Protein Digest (Reduced+Alkylated)**.
- Click **OK** in the **Select Modifications and Profiles** dialog box.
- Mark the **Trypsin** check box under **Enzymes**.

- For this exercise, you use the sequence as is, but you can add modifications and links to sequences as described in *online Help*.
- You can customize the list of available reagents using the Chemical Data Dictionary; see *online Help* for more information.

7 Review parameters on the Mass Matching tab.

- Click the **Mass Matching** tab in the Protein Digest > Match Tolerances section of the Method Editor window.
- Review the parameters.

8 Review the Matching Rules.

- Click the **Matching Rules** tab in the Protein Digest > Match Tolerances section in the Method Editor.
- Mark the **Allow free cysteines (non-reduced condition)** check box.
- Enter 2 for the **Allow missed cleavages up to**.
- Review the other parameters.

9 Save the method for use in Exercise 7.

- Click **Method > Save As**.
- Type the **File name** *iii_NIST_mAb_ProteinDigest.m*, where *iii* represents your initials.
- Click **Save**.

Protein Digest Workflow

Exercise 3. Interactive Protein Digest Sequence Matching

Steps	Detailed Instructions	Comments
10 Start the match search.	<ol style="list-style-type: none">Click Find and Identify > Match Sequences.Select NIST mAb Digest.d.Click Match.	<p><i>Alternate methods:</i></p> <ul style="list-style-type: none">Click  on the Method Editor toolbar.Click Match Sequences on the Method Editor shortcut menu.
11 Review the results.	<ol style="list-style-type: none">Highlight Biomolecule 3 in the Biomolecules window.Click the Biomolecule Identification Results tab which is tabbed with the Biomolecules window.When you open the window, the window displays the results for the first biomolecule that is selected in the Biomolecules window.Select another sequence match result to view by selecting a different biomolecule in the Biomolecules window.	<ul style="list-style-type: none">If the biomolecule was identified, the ID Techniques Applied column contains Sequence Match.
12 View sequence coverage results.	<ol style="list-style-type: none">If necessary, click View > Sequence Coverage Map.Select a different biomolecule in the Biomolecules table to view a different result.	<ul style="list-style-type: none">Amino acids that are matched are either green (MS/MS) or black (MS-only) in a matched sequence.Amino acids that are not matched are gray.A line is added below the AA sequence to display where peptides have been identified.See the online Help for more information.
To view more information.	<p>Click the following items on the Sequence Coverage Map window shortcut menu to view more information about the sequence:</p> <ul style="list-style-type: none">• Applied Modifications• Specified Applied Links• View Digest List	

Protein Digest Workflow

Exercise 3. Interactive Protein Digest Sequence Matching

Steps

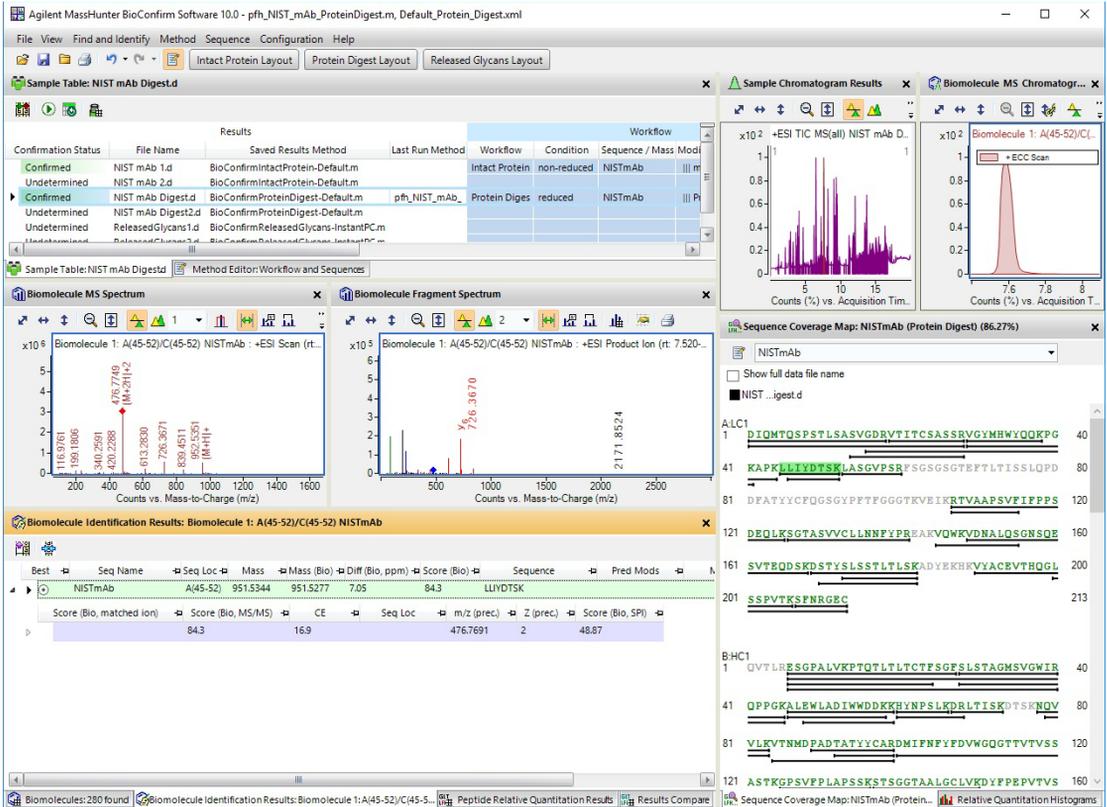
Detailed Instructions

Comments

13 Save the results

a Click **File > Save Results** to save your results to the data file folder.

• You can also click  to save results.



The screenshot displays the Agilent MassHunter BioConfirm Software 10.0 interface. The main window shows a 'Results' table with columns for Confirmation Status, File Name, Saved Results Method, Last Run Method, Workflow, Condition, and Sequence / Mass Mod. The table lists several entries, including 'Confirmed' and 'Undetermined' results for 'NIST mAb Digest.d'. Below the table, there are several panels: 'Biomolecule MS Spectrum' showing a mass spectrum with peaks at m/z 476.2749 and 756.3670; 'Biomolecule Fragment Spectrum' showing a mass spectrum with peaks at m/z 756.3670 and 2171.9524; 'Biomolecule Identification Results' showing a table of results for 'Biomolecule 1: A(45-52)/C(45-52) NISTmAb' with columns for Best, Seq Name, Seq Loc, Mass, Mass (Bio), Diff (Bio, ppm), Score (Bio), Sequence, and Pred Mods; and 'Sequence Coverage Map: NISTmAb (Protein Digest) (86.27%)' showing a protein sequence with coverage bars and a 'Use for %Quant' checkbox checked for the first two rows.

Best	Seq Name	Seq Loc	Mass	Mass (Bio)	Diff (Bio, ppm)	Score (Bio)	Sequence	Pred Mods
1	NISTmAb	A(45-52)	951.5344	951.5277	7.05	84.3	LIVDTSK	
2			84.3					
3			169					
4			476.7891					
5			2					
6			48.87					

14 Review the results in the Peptide Relative Quantitation Results window.

a Click the **Peptide Relative Quantitation Results** tab. It is tabbed with the Biomolecules window.
 b Click the first triangle next to the first row.
 c Note that **Use for %Quant** is marked for both rows.

• For Protein Digest workflow, the software uses the **Use for %Quant** check box in the Peptide Relative Quantitation Results window.

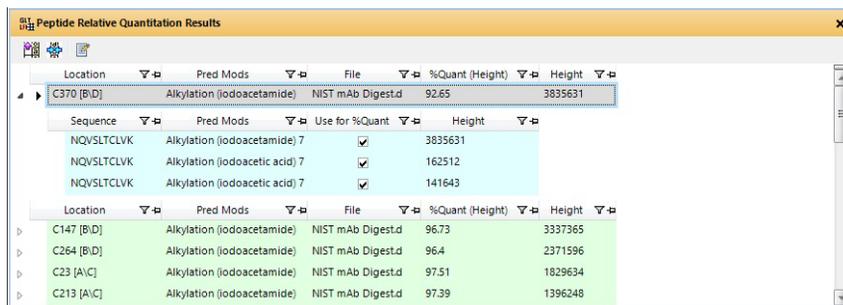
Protein Digest Workflow

Exercise 3. Interactive Protein Digest Sequence Matching

Steps

Detailed Instructions

Comments

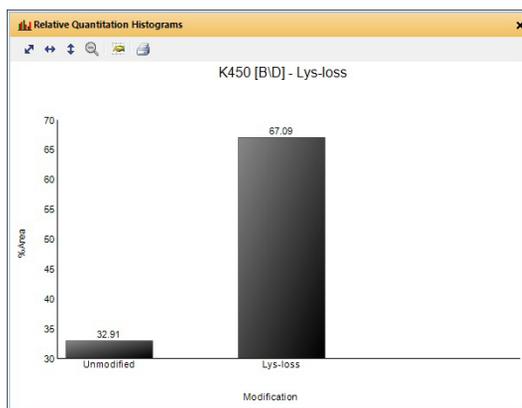


Location	Pred Mods	File	%Quant (Height)	Height
C370 [B:D]	Alkylation (iodoacetamide)	NIST mAb Digest.d	92.65	3835631
Sequence				
NQVSLTCLVK	Alkylation (iodoacetamide) 7		<input checked="" type="checkbox"/>	3835631
NQVSLTCLVK	Alkylation (iodoacetic acid) 7		<input checked="" type="checkbox"/>	162512
NQVSLTCLVK	Alkylation (iodoacetic acid) 7		<input checked="" type="checkbox"/>	141643
Location				
C147 [B:D]	Alkylation (iodoacetamide)	NIST mAb Digest.d	96.73	3337365
C264 [B:D]	Alkylation (iodoacetamide)	NIST mAb Digest.d	96.4	2371596
C23 [A:C]	Alkylation (iodoacetamide)	NIST mAb Digest.d	97.51	1829634
C213 [A:C]	Alkylation (iodoacetamide)	NIST mAb Digest.d	97.39	1396248

15 Review the results in the Relative Quantitation Histograms window.

- Click the **Relative Quantitation Histograms** window.
- Click the first triangle next to the first row.
- Note that **Use for %Quant** is marked for both rows.

- For Protein Digest workflow, the software uses the **Use for %Quant** check box in the Peptide Relative Quantitation Results window.



16 Repeat the interactive processing with *NIST mAb Digest2.d*.

- Open the data file **NIST mAb Digest2.d** (see **step 3**).
- Select Find Peptides in the Method Editor and verify the parameters (**step 4**).
- Find biomolecules (**step 5**).
- Match sequences (**step 10**).
- Save the results to the NIST mAb Digest2.d file (**step 13**).

- Most of the processing parameters used for the first data file are the same for the second data file.
- These results are used in **“Exercise 8. Using Result Review mode”** on page 36.

Exercise 4. Automated Protein Digest Workflow

This exercise guides you through the setup of a worklist to automatically confirm the presence of NIST mAb in a previously acquired sample.

If you select the Protein Digest workflow, the Find Peptides algorithm runs and uses the enzymes selected in the Workflow and Peptides section and then runs the Protein Digest matching rules.

Steps	Detailed Instructions	Comments
1 Open the method.	<ol style="list-style-type: none"> a Click Method > Open. b Select the <i>iii_NIST_mAb_ProteinDigest.m</i> folder. c Click Open. 	<ul style="list-style-type: none"> • This method was created in Exercise 3 (<i>iii</i> represents your initials).
2 Display the Method Automation > Workflow and Sequences section in the Method Editor.	<ol style="list-style-type: none"> a If the Method Editor is not visible, click View > Method Editor. b Click Method Automation > Workflow and Sequences in the Method Editor window. 	<ul style="list-style-type: none"> • You can instead click the Method Editor button, , on the main toolbar.
3 Select the appropriate workflow.	<ol style="list-style-type: none"> a Select Protein Digest for the Workflow. b Select the Condition. c Verify that NISTmAb is the sequence. d Verify that Protein Digest (Reduced+Alkylated) is the Mods and Profiles. e Mark the Trypsin check box. 	<ul style="list-style-type: none"> • The Protein Digest workflow automatically runs the following actions: <ul style="list-style-type: none"> • Find Peptides • Match Sequences
4 Save the method.	<ul style="list-style-type: none"> • Click Method > Save. 	
5 Run the method workflow or run method automation.	<ul style="list-style-type: none"> • Click Method > Run Method Workflow. • Click Method > Run Method Automation (Workflow + Reports). • Click  on the Method Editor toolbar. • Reprocess the sample. See “Exercise 7. Reprocessing Samples” on page 33. 	<ul style="list-style-type: none"> • Method Automation first runs the method workflow and then extracts additional chromatograms, generates a biomolecule report, and exports results.
6 (optional) Review the printed Biomolecule reports.	<ul style="list-style-type: none"> • If you clicked Run Method Automation (Workflow + Reports), then a report is generated automatically. • You can click File > Print > Biomolecule Report to generate a report for the current sample. 	<ul style="list-style-type: none"> • You set report options in the Method Editor window in the Method Automation > Reports section. • If you are creating a report interactively, you specify some of these parameters in the Print Biomolecule Report dialog box.

Released Glycans Workflow

Exercise 4. Automated Protein Digest Workflow

Released Glycans Workflow

The steps outlined below show the workflow for Released Glycans with Agilent MassHunter BioConfirm Software.

Step 1 - Open the data file of interest.

Step 2 - Open a BioConfirm Method or create a new one.

Step 3-Select the **Workflow** on the **Workflow and Sequences** tab.

Step 4 - Select the **Target glycan source**.

Step 5 - Select the tag which you used. 2-AB and InstantPC are listed, and you can create your own.

Step 6 - Run the Method Workflow.

Step 9 - Review the results which are shown in these windows:

Sample Chromatogram Results

Biomolecule MS Chromatogram

Biomolecules table

Biomolecule Identification Results

Biomolecule MS Spectrum

Biomolecule Fragment Spectrum

Glycan Structure Viewer

Results Compare

Relative Quantitation Histograms

Step 9 - Print report.

Exercise 5. Interactive Released Glycans

This exercise shows you how to find released glycans interactively.

If you select the Released Glycans workflow, the Find Glycans algorithm runs. See “**Before you start**” on page 2.

Steps	Detailed Instructions	Comments
1 Open the method to use as a starting point for the new method.	<ol style="list-style-type: none"> a Click Method > Open. b Select the BioConfirmReleasedGlycans-Default.m folder. c Click Open. 	<ul style="list-style-type: none"> • The parameters in the BioConfirmProteinDigest-Default.m method are a good starting point for Protein Digests.
2 Open the example sample file.	<ol style="list-style-type: none"> a Click File > Open Data File. b Locate the ReleasedGlycans1.d folder. c Click Open. 	<ul style="list-style-type: none"> • The TIC is automatically displayed in the Sample Chromatogram Results window.
3 Load the Released Glycans layout.	<ul style="list-style-type: none"> • Click Released Glycans Layout in the main toolbar. 	
4 Review the parameters in the Find Peptides section in the Method Editor window.	<ol style="list-style-type: none"> a Select Method Automation > Workflow and Sequences in the Method Editor window. b Enter <code>Example</code> in the Glycan group. c Clear Require RT match if database contains a RT for the target glycan. d Select Released Glycans > Find Glycans in the Method Editor window. e Select <code>Glycans_mAb_AM_PCD.cdb</code> for the Target glycan source. f Click the Tag tab. g Click the option for the correct tag. For the example data file, click InstantPC. 	<ul style="list-style-type: none"> • You can change the default parameters as described in the next steps. • For some data files, you will need to use different parameters as described in the <i>online Help</i>. • A very low peak height filter can result in greater sequence coverage but requires much more time to process.
5 Find biomolecules.	<ol style="list-style-type: none"> a Click  on the Method Editor toolbar to start the biomolecule search. b When processing is complete, review the results in the Biomolecules window. c Click View > Glycan Structure Viewer. 	
6 Save the method for use in Exercise 6.	<ol style="list-style-type: none"> a Click Method > Save As. b Type the File name <code>iii_ReleasedGlycans_InstantPC.m</code>, where <i>iii</i> represents your initials. c Click Save. 	

Released Glycans Workflow

Exercise 5. Interactive Released Glycans

Steps

7 Review the results.

Detailed Instructions

- In the Biomolecules window, click the header of the **Area (Glycan)** column to sort the table by this column. If necessary, click the header again so that the largest areas are at the top of the table.
- Highlight **G0F** in the Biomolecules window.
- Click the **Biomolecule Identification Results** tab. The Biomolecules Identification Results window is tabbed with the Biomolecules window.
- When you open the window, the window displays the results for the first biomolecule that is selected in the Biomolecules window.

Comments

- Several changes were made to the default layout for the image below. The Glycan Structure Viewer window is visible. Also, the **Flags (Tgt)** column was moved.
- The Relative Quantitation Histograms window only contains information when you run a workflow.

The screenshot displays the Agilent MassHunter BioConfirm Software 10.0 interface. The main window is titled 'Released Glycans' and shows the following components:

- Results Table:** A table with columns for Confirmation Status, File Name, Saved Results Method, Workflow, Condition, and Sequence. The 'G0F' entry is highlighted.
- Sample Chromatogram Results:** A plot showing the Total Ion Chromatogram (TIC) scan for the released glycans.
- Biomolecule MS Chromatogram:** A plot showing the mass spectrum for the selected biomolecule (G0F) at 14.98 minutes.
- Biomolecule MS Spectrum:** A plot showing the mass spectrum for the selected biomolecule (G0F) at 14.98 minutes.
- Glycan Structure Viewer: G0F:** A diagram showing the chemical structure of the G0F glycan.
- Biomolecules: 145 found:** A table listing 145 biomolecules with columns for Name, Notes, Flags (Tgt), Diff (Tgt, ppm), Score (Tgt), Hits, Use for %Quant, and Area (Gly). The 'G0F' entry is highlighted.
- Relative Quantitation Histograms:** A window showing no data to display.

8 Save the results

- Click **File > Save Results** to save your results to the data file folder.

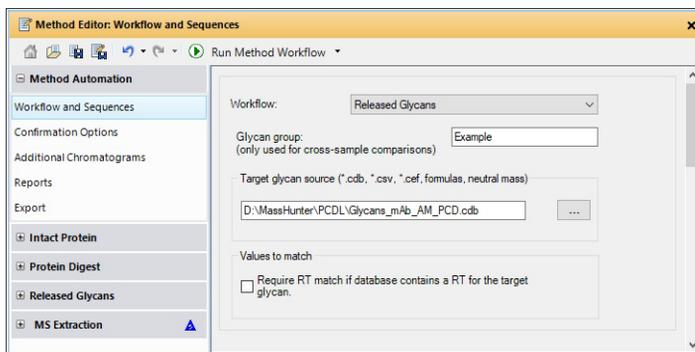
- You can also click  to save results.

Exercise 6. Automated Released Glycans Workflow

This exercise guides you through the setup of a worklist to automatically run the Released Glycans workflow.

If you select the Released Glycans workflow, the Find Glycans algorithm runs and uses the **Target glycan source** selected in the Workflow and Peptides section.

Steps	Detailed Instructions	Comments
1 Open the method.	<ol style="list-style-type: none"> Click Method > Open. Select the <i>iii_ReleasedGlycans_InstantPC.m</i> folder. Click Open. 	<ul style="list-style-type: none"> This method was created in Exercise 5 (<i>iii</i> represents your initials).
2 Display the Method Automation > Workflow and Sequences section in the Method Editor.	<ol style="list-style-type: none"> If the Method Editor is not visible, click View > Method Editor. Click Method Automation > Workflow and Sequences in the Method Editor window. 	<ul style="list-style-type: none"> You can instead click the Method Editor button, , on the main toolbar.
3 Select the appropriate workflow.	<ol style="list-style-type: none"> Select Released Glycans for the Workflow. Enter <i>Example</i> in the Glycan group. Select Glycans_mAb_AM_PCD.cdb as the Target glycan source. Clear the Require RT match if database contains a RT for the target glycan check box. 	<ul style="list-style-type: none"> The Released Glycans workflow automatically runs the Find Glycans algorithm. The Glycan group is used to organize the results in the Results Compare window.
4 Save the method.	<ul style="list-style-type: none"> Click Method > Save. 	



Released Glycans Workflow

Exercise 6. Automated Released Glycans Workflow

Steps	Detailed Instructions	Comments
5 Run the method workflow or run method automation.	<ul style="list-style-type: none">Click Method > Run Method Workflow.Click Method > Run Method Automation (Workflow + Reports).Click  on the Method Editor toolbar when the Workflow and Sequences section is showing.Reprocess the sample. See “Exercise 7. Reprocessing Samples” on page 33.	<ul style="list-style-type: none">Method Automation first runs the method workflow and then extracts additional chromatograms, generates a biomolecule report, and exports results.The Workflow column is set to “Released Glycans”. The Relative Quantitation Histograms window and the Results Compare window contain results.
6 (optional) Review the printed Biomolecule reports.	<ul style="list-style-type: none">If you clicked Run Method Automation (Workflow + Reports), then a report is generated automatically.You can click File > Print > Biomolecule Report to generate a report for the current sample.	<ul style="list-style-type: none">You set report options in the Method Editor window in the Method Automation > Reports section.If you are creating a report interactively, you specify some of these parameters in the Print Biomolecule Report dialog box.

Released Glycans Workflow

Exercise 6. Automated Released Glycans Workflow

Steps

Detailed Instructions

Comments

The screenshot displays the Agilent MassHunter BioConfirm Software interface for the Released Glycans Workflow. The main window is titled "Agilent MassHunter BioConfirm Software 10.0 - pfh_ReleasedGlycans_InstanPC.m, Default_Released_Glycans.xml".

Results Table:

Confirmation Status	File Name	Saved Results Method	Workflow	Condition	Sequence / N
Confirmed	NIST mAb 1.d	BioConfirmIntactProtein-Default.m	Intact Protein	non-reduced	NISTmAb
Undetermined	NIST mAb 2.d	BioConfirmIntactProtein-Default.m	Intact Protein	non-reduced	NISTmAb
Confirmed	NIST mAb Digest.d	BioConfirmProteinDigest-Default.m	Protein Digest	reduced	NISTmAb
Confirmed	NIST mAb Digest2.d	BioConfirmProteinDigest-Default.m	Protein Digest	reduced	NISTmAb
Undetermined	ReleasedGlycans1.d	pfh_ReleasedGlycans_InstanPC.m	Released Glyc		
Undetermined	ReleasedGlycans2.d	BioConfirmReleasedGlycans-InstanPC.m	Released Glyc		

Sample Chromatogram Results: Shows a Total Ion Chromatogram (TIC) scan for the released glycan sample, with a peak at 14.468 minutes.

Biomolecule MS Chromatogram: Shows a chromatogram for the biomolecule GOF (C70 H113 N7 O42) with a peak at 14.98 minutes.

Biomolecule MS Spectrum: Shows the mass spectrum for the biomolecule GOF, with a base peak at m/z 1497.

Glycan Structure Viewer: Displays the chemical structure of the glycan GOF, which is a branched N-glycan with a core fucose.

Biomolecules: 145 found:

Name	Notes	Flags (Tgt)	Diff (Tgt, ppm)	Score (Tgt)	Hits	Use for %Quant	Area (Gly)
GOF	Hex3HexNAc4dH	multiple IDs	1.11	99.1	2	<input checked="" type="checkbox"/>	2910586
G1F	Hex4HexNAc4dH	multiple IDs	0.7	99.02	2	<input checked="" type="checkbox"/>	1236757
G1F	Hex4HexNAc4dH	multiple IDs	0.84	98.49	2	<input checked="" type="checkbox"/>	5980375
H4N4S1	Hex4HexNAc4Ne		-5.06	82.48	1	<input checked="" type="checkbox"/>	1645265
HSN4F1	Hex5HexNAc4dH		0.21	99.81	1	<input checked="" type="checkbox"/>	1315894
GOF-GlcNAc	Hex3HexNAc3dH	multiple IDs	0.51	99.42	2	<input checked="" type="checkbox"/>	1302664
HSNSF1S2	Hex5HexNAc5dH	low score	7.93	57.94	1	<input checked="" type="checkbox"/>	7164260
H4N4S1	Hex4HexNAc4Ne		-5.39	81.23	1	<input checked="" type="checkbox"/>	6330075
HSN3S1	Hex5HexNAc3He		-0.75	98.36	2	<input checked="" type="checkbox"/>	5181136
H4N3S1	Hex4HexNAc3He		-5.86	79.74	1	<input checked="" type="checkbox"/>	4966014
H3NSF1	Hex3HexNAc5dH		-0.04	99.86	1	<input checked="" type="checkbox"/>	4178319
HSN4F1	Hex5HexNAc4dH		-0.08	99.79	1	<input checked="" type="checkbox"/>	4097144

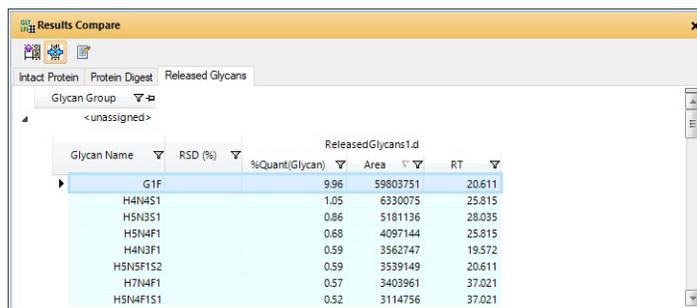
Results Compare:

Glycan Name	RSD (%)	%Quant(Glycan)	Area	RT
H6NF1S2	0	0	9090	51.19
H6NF1S1	0	0	12858	50.93
H6NF5S1	0.04	0.04	251939	38.1
H6NF4F1S1	0.01	0.01	69946	37.30
HSN4F2S1	0	0	2996	40.66
HSN5S1	0.03	0.03	152807	34.77
HSN4F1S1	0.52	0.52	3114756	37.02
HSN3S1	0.86	0.86	5181136	28.03
H6NSF2S3	0.04	0.04	226037	50.75

Released Glycans Workflow

Exercise 6. Automated Released Glycans Workflow

Steps	Detailed Instructions	Comments
7 Review the results in the Results Compare window.	<ol style="list-style-type: none">Click the Results Compare window.Click the Released Glycans tab.Note that RSD (%) is empty because only one sample is selected. If you select two or more samples and they belong to the same Glycan Group, then the results are shown in the same table. If the same glycan is in multiple samples, then the RSD (%) is calculated.	<ul style="list-style-type: none">For Released Glycans workflow, the software uses the Use for %Quant check box in the Biomolecules window.



Glycan Name	RSD (%)	%Quant(Glycan)	Area	RT
G1F		9.96	59803751	20.611
H4N4S1		1.05	6330075	25.815
H5N3S1		0.86	5181136	28.035
H5N4F1		0.68	4097144	25.815
H4N3F1		0.59	3562747	19.572
H5N5F1S2		0.59	3539149	20.611
H7N4F1		0.57	3403961	37.021
H5N4F1S1		0.52	3114756	37.021

8 Save the results and close the data file.	<ol style="list-style-type: none">Click File > Save Results.Click File > Close All.Click Yes to save results.
---	---

Review Results

Exercise 7. Reprocessing Samples

This exercise shows you how to reprocess samples in the Sample Table. You can quickly check the Confirmation Status of each sample and determine if you need to reprocess the sample.

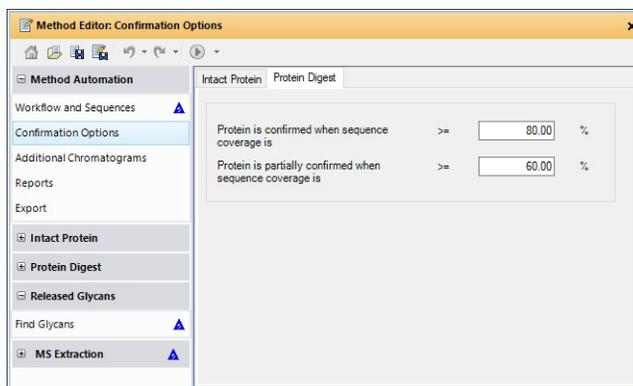
Steps	Detailed Instructions	Comments
1 Open several data files.	<p>a Click File > Open Sample Files.</p> <p>b Select these example files: NIST mAb1.d, NIST mAb 2.d, NIST mAb Digest.d, NIST mAb Digest2.d, ReleasedGlycans1.d, and ReleasedGlycans2.d.</p> <p>c Mark the Load result data check box.</p> <p>d Click Open.</p>	<ul style="list-style-type: none"> To select multiple files, click the first file. Then, press Shift and click the last file.
2 Review results in the Sample Table window.	<p>a Look at the Confirmation Status column.</p>	<ul style="list-style-type: none"> If you saved results, the table contains information on confirmation.

Results		Workflow					
Confirmation Status	File Name	Workflow	Condition	Sequence / Mass	Modification	Enzyme	Glycan Group
Confirmed	NIST mAb 1.d	Intact Protein	non-reduced	NISTmAb	mAb		
Undetermined	NIST mAb 2.d	Intact Protein	reduced				
Confirmed	NIST mAb Digest.d	Protein Digest	reduced	NISTmAb	Protein Digest (Reduced+Alkylated)	Trypsin,LysC	
Confirmed	NIST mAb Digest2.d	Protein Digest	reduced	NISTmAb	Protein Digest (Reduced+Alkylated)	Trypsin,LysC	
Undetermined	ReleasedGlycans1.d	Released Glycans					Example
Undetermined	ReleasedGlycans2.d	Intact Protein	reduced				

Review Results

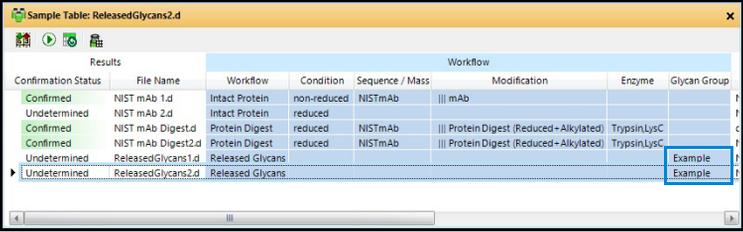
Exercise 7. Reprocessing Samples

Steps	Detailed Instructions	Comments
3 Review values in Method Automation > Confirmation Options.	<ol style="list-style-type: none">Click View > Method Editor, if necessary.Select Method Automation > Confirmation Options.Click the Intact Protein tab.Review selection for the Intact Protein match found but not for the most abundant peak option.Click the Protein Digest tab.Review selection for the Protein is partially confirmed when sequence coverage is >= option.	<ul style="list-style-type: none">These tabs explain what it means to be Confirmed and Partially confirmed.You are not changing these options. You are only seeing what the software checks to determine if the protein is confirmed.



Review Results

Exercise 7. Reprocessing Samples

Steps	Detailed Instructions	Comments
4	<p>Reprocess the ReleasedGlycans2.d data file.</p> <ol style="list-style-type: none"> In the Sample Table, click the row containing ReleasedGlycans2.d. Click Method > Open. Select the <i>iii_ReleasedGlycans-InstantPC.m</i> folder. Click Open. Click the  button to open the Reprocess Sample dialog box. Select Released Glycans for the workflow. Enter Example for the Glycan group. Select Glycans_mAb_AM_PCD.cdb for the Target glycan source. Click Reprocess. 	<ul style="list-style-type: none"> To reprocess a sample, you need to first load the correct method and then complete the Reprocess Sample dialog box. You can also double-click the Sample Table row to open the Reprocess Sample dialog box. You can either use the current method, or if you have previously saved results, you can use the sample result method.
	<p>Both data files that used the Released Glycans workflow have the same Glycan Group.</p>	
		
5	<p>Save the results for the samples that you reprocessed.</p> <ol style="list-style-type: none"> Click File > Save Results. Click Save. 	

Review Results

Exercise 8. Using Result Review mode

Exercise 8. Using Result Review mode

This exercise shows you how to use the Result Review mode. When this mode is enabled, you cannot edit a method. You also cannot run the algorithms in the Find and Identify menu.

Steps	Detailed Instructions	Comments
1 Enable Result Review mode.	<ul style="list-style-type: none">Click Configuration > Enable Result Review (Disables Method Editing).	<ul style="list-style-type: none">You can toggle this mode off by clicking this same comment again.
2 Review results in Sample Table window.	<ul style="list-style-type: none">All of the options in this window are available except for the Run Method Workflow button. You can still reprocess samples.	

Confirmation Status	Results		Workflow					
	File Name	Saved Results Method	Workflow	Condition	Sequence / Mass	Modification	Enzyme	Glycan Group
Confirmed	NIST mAb 1.d	pft_ReleasedGlycans_InstanPC.m	Intact Protein	non-reduced	NISTmAb	mAb		
Undetermined	NIST mAb 2.d	pft_ReleasedGlycans_InstanPC.m	Intact Protein	reduced				
Confirmed	NIST mAb Digest.d	pft_ReleasedGlycans_InstanPC.m	Protein Digest	reduced	NISTmAb	Protein Digest (Reduced+Alkylated)	Trypsin,LysC	
Confirmed	NIST mAb Digest2.d	pft_ReleasedGlycans_InstanPC.m	Protein Digest	reduced	NISTmAb	Protein Digest (Reduced+Alkylated)	Trypsin,LysC	
Undetermined	ReleasedGlycans1.d	pft_ReleasedGlycans_InstanPC.m	Released Glycans					Example
Undetermined	ReleasedGlycans2.d	pft_ReleasedGlycans_InstanPC.m	Released Glycans					Example

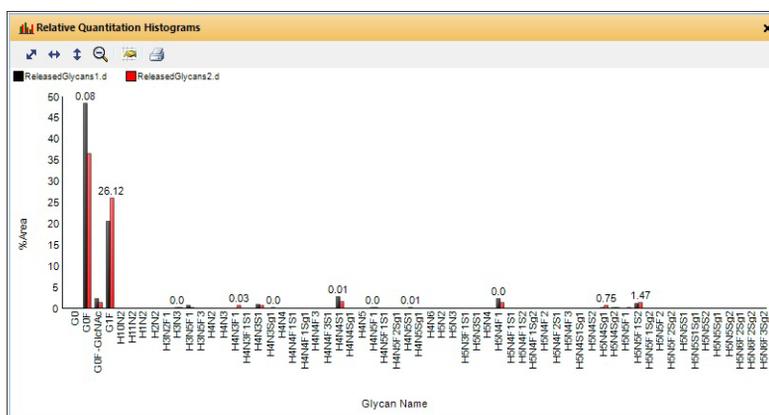
Review Results

Exercise 8. Using Result Review mode

Steps	Detailed Instructions	Comments
3	<p>Review and compare the <i>ReleasedGlycans1.d</i> and <i>ReleasedGlycans2.d</i> data files.</p> <p>a In the Sample Table, click the row containing ReleasedGlycans1.d.</p> <p>b Press the Shift key and click the row containing ReleasedGlycans2.d.</p> <p>c Review the results in the Results Compare window. The RSD (%) column has been calculated for glycans that are in both samples.</p>	

Glycan Name	RSD (%)	ReleasedGlycans1.d			ReleasedGlycans2.d		
		%Quant(Glycan)	Area	RT	%Quant(Glycan)	Area	RT
G1F	9.5	9.96	59803751	20.611	9.24	52275585	20.602
H4N4S1	105	1.05	6330075	25.815	7.58	42841859	25.807
H5N3S1	1.4	0.86	5181136	28.035	0.9	5080633	28.018
H5N4F1	77	0.68	4097144	25.815	0.21	1209200	26.58
H4N3F1	59.5	0.59	3562747	19.572	0.26	1453310	20.595
H5N5F1S2	57.1	0.59	3539149	20.611	1.47	8335163	19.564
H7N4F1	4.9	0.57	3403961	37.021	0.56	3174200	37.021
H5N4F1S1	0.6	0.52	3114756	37.021	0.55	3086343	37.021
H5N2	7.1	0.49	2959099	16.438	0.47	2677458	16.447
H4N5F1	35	0.42	2523553	22.723	0.27	1523201	22.715
G0F-GlcNAc	93.6	0.32	1931147	19.572	1.68	9900305	14.992
H5N4S1Sg1	9.7	0.3	1807009	39.465	0.28	1576053	39.473
H4N4F1Sg1	43.6	0.26	1586348	30.662	0.15	839075	30.645
H5N4S2	3.9	0.23	1356469	39.473	0.23	1282857	39.482

- d Review the Relative Quantitation Histograms window.
- You can visually compare the relative quantitation results for different glycans.



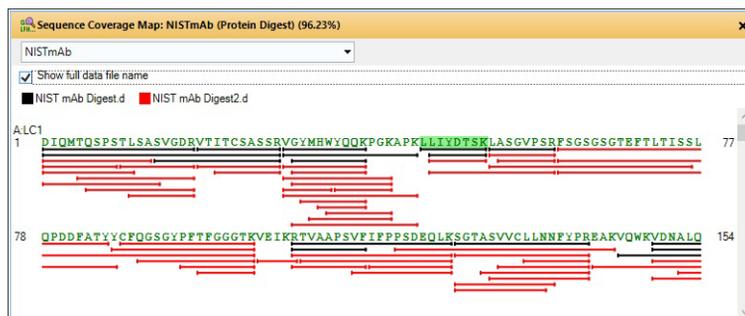
Review Results

Exercise 8. Using Result Review mode

Steps	Detailed Instructions	Comments
4	<p>Review and compare the <i>NIST mAb Digest.d</i> and <i>NIST mAb Digest2.d</i> data files.</p> <ol style="list-style-type: none"> In the Sample Table, click the row containing NIST mAb Digest.d. Press the Shift key and click the row containing NIST mAbDigest2.d. Click the Protein Digest Layout button in the main toolbar. Review the results in the Results Compare window. The RSD (%) column has been calculated for many modifications that are in both samples. 	•

NISTmAb				NIST mAb Digest.d		NIST mAb Digest2.d	
Location	Modification	Average (Height)	%RSD (Height)	%Quant (Height)	Height	%Quant (Height)	Height
C370 [B/D]	Alkylation (Iodo...)	2124028	114	92.65	3835631	73.3	412425
C147 [B/D]	Alkylation (Iodo...)	1895628	107.6	96.73	3337365	91.07	453891
C264 [B/D]	Alkylation (Iodo...)	1204864	136.9	96.4	2371596	58.96	38132
C23 [A/C]	Alkylation (Iodo...)	1085689	96.9	97.51	1829634	100	341744
C213 [A/C]	Alkylation (Iodo...)	726794	130.3	97.39	1396248	100	57340
C97 [B/D]	Alkylation (Iodo...)	306169	125.4	100	577691	100	34648
C193 [A/C]	Alkylation (Iodo...)	423883	35.3	61.9	318141	100	529624
M255 [B/D]	Oxidation (M)	198278	52.9	12.41	272404	15	124153
C428 [B/D]	Alkylation (Iodo...)	368817	18.9	74.04	319406	87.48	418228
C370 [B/D]	Alkylation (Iodo...)	224645	60.1	7.35	304155	25.70	145134

- Review the results in the Sequence Coverage Map window. A legend is added to the top of the window and the lines under the sequence are color coded to show which sample is described.



Review Results

Exercise 8. Using Result Review mode

Steps	Detailed Instructions	Comments
	<p>f Review the results in the Peptide Relative Quantitation Results window. The Location C147 [B/D] has two rows because this location is in both samples.</p>	<ul style="list-style-type: none"> For Protein Digest workflows, you mark the Use for %Quant check box in the Peptide Relative Quantitation Results window.

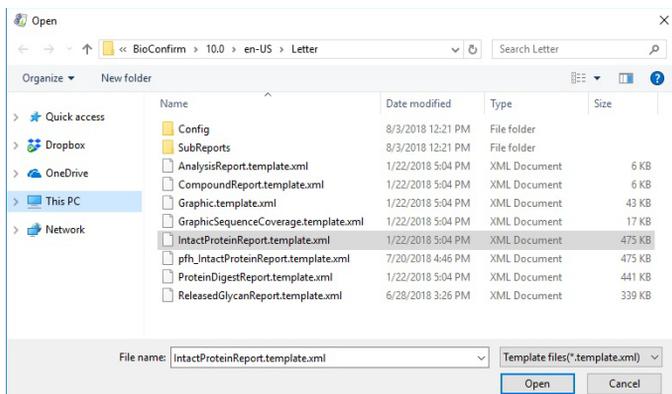
Location	Pred Mods	File	%Quant (Height)	Height
C133 [A/C]	Alkylation (iodoacetic acid)	NIST mAb Digest.d	100	99284
C133 [A/C]	Alkylation (iodoacetamide)	NIST mAb Digest2.d	73.29	136442
C133 [A/C]	Alkylation (iodoacetic acid)	NIST mAb Digest2.d	26.71	49714
C147 [B/D]	Alkylation (iodoacetic acid)	NIST mAb Digest.d	96.73	3337365
C147 [B/D]	Alkylation (iodoacetic acid)	NIST mAb Digest.d	3.27	112950
Sequence				
STSGGTAALGCLVK	Alkylation (iodoacetamide) 1	<input checked="" type="checkbox"/>	3337365	
LGCLVK	Alkylation (iodoacetamide) 3	<input type="checkbox"/>	187186	
STSGGTAALGCLVK	Alkylation (iodoacetic acid) 1	<input checked="" type="checkbox"/>	112950	
SGGTAALGCLVK	Alkylation (iodoacetamide) 9	<input type="checkbox"/>	26054	
Location				
C147 [B/D]	Alkylation (iodoacetamide)	NIST mAb Digest2.d	91.07	453891
Sequence				
STSGGTAALGCLVK	Alkylation (iodoacetamide) 1	<input checked="" type="checkbox"/>	278669	
LGCLVK	Alkylation (iodoacetamide) 3	<input checked="" type="checkbox"/>	158833	
STSGGTAALGCLVK	Alkylation (iodoacetic acid) 1	<input checked="" type="checkbox"/>	28345	
GGTAALGCLVK	Alkylation (iodoacetic acid) 8	<input checked="" type="checkbox"/>	12788	
AALGCLVK	Alkylation (iodoacetic acid) 5	<input checked="" type="checkbox"/>	12047	
SGGTAALGCLVK	Alkylation (iodoacetic acid) 9	<input checked="" type="checkbox"/>	4343	
STSGGTAALGCLVK		<input checked="" type="checkbox"/>	3378	
Location				
C147 [B/D]	Alkylation (iodoacetic acid)	NIST mAb Digest2.d	8.25	41133
C193 [A/C]	Alkylation (iodoacetamide)	NIST mAb Digest.d	61.9	318141
C193 [A/C]	Alkylation (iodoacetic acid)	NIST mAb Digest.d	38.1	195816
C193 [A/C]	Alkylation (iodoacetamide)	NIST mAb Digest2.d	100	529624

- 5 Save the results for the samples that you reprocessed.
- a Click **File > Save Results**.
- b Click **Save**.

Exercise 9. Using Report Builder

This exercise shows you the program to modify PDF templates. If you click **Use PDF Report Builder** in the Method Automation > Reports > Templates tab, then you can use Report Builder to modify those templates.

Steps	Detailed Instructions	Comments
1 Open Report Builder program.	<ol style="list-style-type: none">Double-click Report Builder in the Tools for BioConfirm 10.0 folder in the Agilent MassHunter Workstation program folder.In Windows 10, click Agilent MassHunter Report Builder > Report Builder 10.0.	<ul style="list-style-type: none">You can also start the Report Builder program when you click one of the Edit button next to a template in the Method Automation > Reports > Template tab. These Edit buttons are only available if you click Configuration > Show Advanced Settings.
2 Open an existing template.	<ol style="list-style-type: none">Click File > Open > Browse.Select a template and click Open.	<ul style="list-style-type: none">Report templates are installed in the D:\MassHunter\Report Templates\BioConfirm folder.

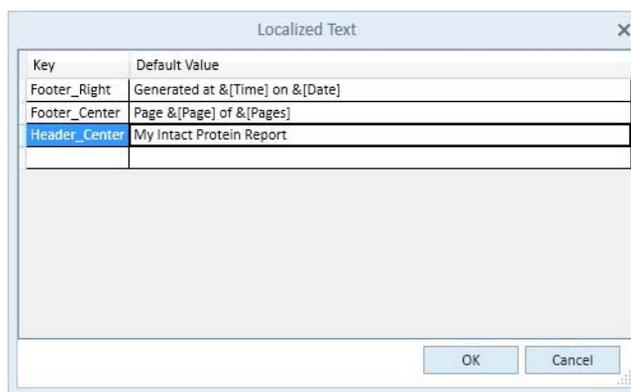


Agilent recommends that you do not modify the default templates. Instead, make a copy of the template and modify the copy.

Review Results

Exercise 9. Using Report Builder

Steps	Detailed Instructions	Comments
3	Review the template in Report Builder. a Click an item in the template. Notice that the right pane changes. b Click the title of the report. c In the right pane, click Localized Text in the Content section. d Click the ... button. The Localized Text dialog box opens. e Click the Header_Center . f Enter My Intact Protein Report . g Click OK .	<ul style="list-style-type: none">• The left pane shows the template. The right pane shows the parameters for the current selection.• You can make many different changes to the report. This exercise only shows you one possibility. Press F1 to access the online Help to learn more about customizing a report template.



Review Results

Exercise 9. Using Report Builder

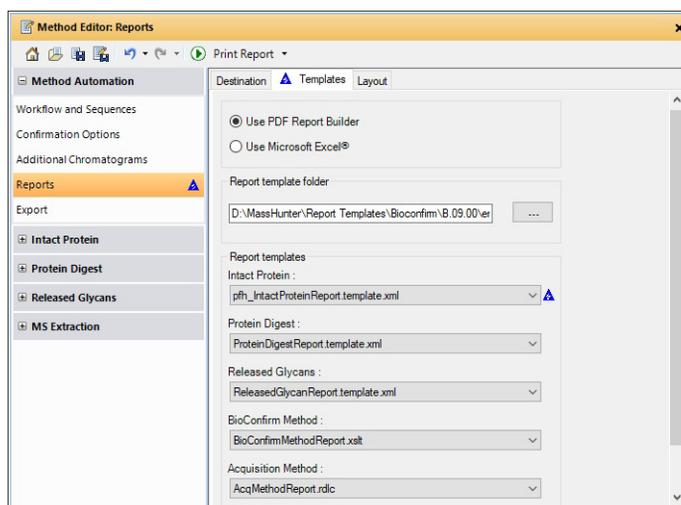
Steps	Detailed Instructions	Comments
4 Save the template.	<ol style="list-style-type: none"> Click File > Save As > Browse. Enter a file name and click Save. Close the Report Builder program. 	<ul style="list-style-type: none"> You can instead click File > Save, and the file is saved to the current method. Agilent recommends that you do not modify the default templates.

The screenshot shows the Agilent Report Builder software interface. The window title is "pfh_IntactProteinReport.template - Report Builder (BioConfirm 10.0 Intact Protein Report)". The interface includes a ribbon with "File", "Home", "Edit", "Font", "Alignment", "Colors", "Page", "Textbox", "Image", "List", "Table", and "Insert" tabs. The main workspace displays a report template for "{L:My Intact Protein Report}" with sections for "GlobalReportProperties", "Sample Information", "Additional Walkup Information", "Matched Sequences", and "Sample Chromatogram List". A right-hand pane shows "Page1" settings, including "General" (ID, Page, Page Orientation, Type) and "Margin" (Margins: 0.75in, 0.35in, 0.25in, 0.25in).

Review Results

Exercise 9. Using Report Builder

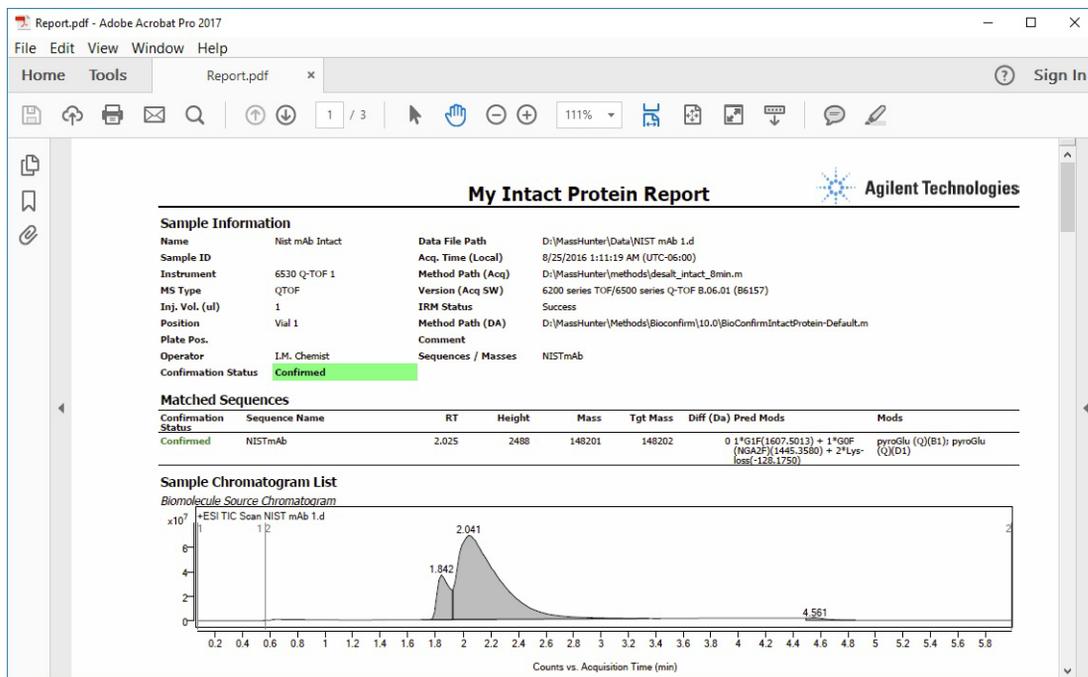
Steps	Detailed Instructions	Comments
5	Use this new template in a method. <ol style="list-style-type: none">Open the Method Editor window. Click View > Method Editor if it is not visible.Select Method Automation > Reports.Click the Templates tab.Select the changed report for the corresponding report template type. In this example, the Intact Protein report template was modified.	<ul style="list-style-type: none">Different reports use different report templates. If you modified an Intact Protein report template, then you select the modified template for the Intact Protein report template.When you print a biomolecule report, the report template corresponding to the selected workflow is used.



Review Results

Exercise 9. Using Report Builder

Steps	Detailed Instructions	Comments
6 Print a Biomolecule report.	<ol style="list-style-type: none"> Click File > Print > Biomolecule Report. Mark the Print preview check box. Click OK. 	<ul style="list-style-type: none"> When you print a biomolecule report, the report template corresponding to the selected workflow is used.



Deconvolution

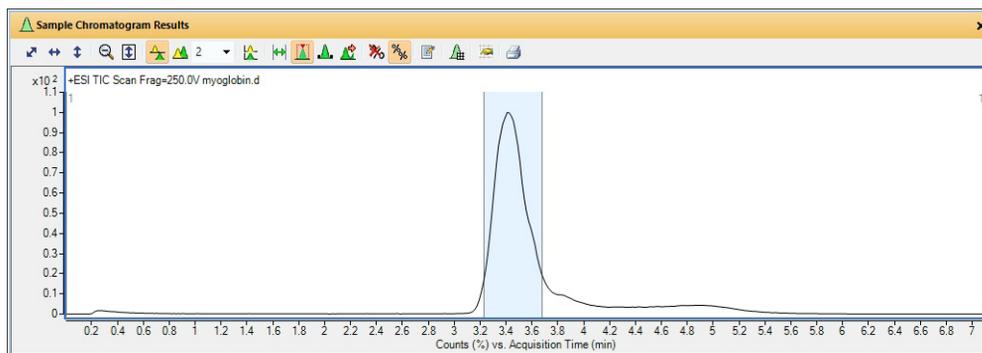
Exercise 10. Interactive Protein Molecular Weight Determination

Deconvolution

Exercise 10. Interactive Protein Molecular Weight Determination

This exercise shows you how to open a data file, extract spectra, deconvolute and view results. Deconvolution software does charge state deconvolution of mass spectra of large molecules with high charge states, such as proteins. See “**Before you start**” on page 2.

Steps	Detailed Instructions	Comments
1 Open the data file.	<ol style="list-style-type: none">Click File > Open Data File.Locate the myoglobin.d folder.Clear the Load Result check box.Click Open.	<ul style="list-style-type: none">The TIC is automatically displayed in the Sample Chromatogram Results window.
2 Extract a peak spectrum.	<ol style="list-style-type: none">Select a range around the peak at 3.5 minutes.Double-click this range.	<ul style="list-style-type: none">To select a range, click one side of the peak and drag to the other side of the peak.



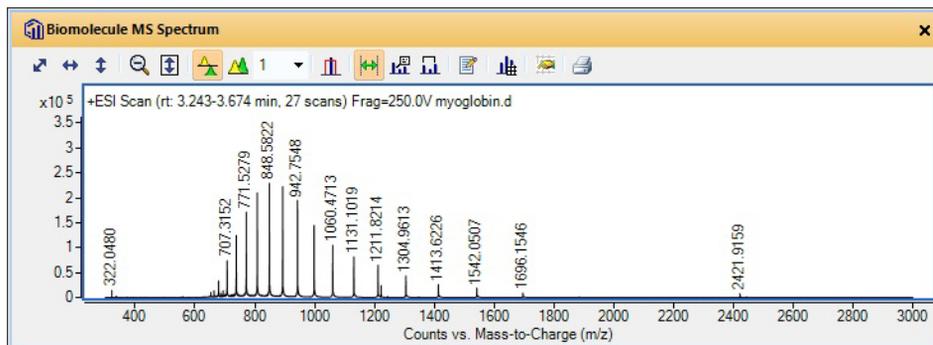
Deconvolution

Exercise 10. Interactive Protein Molecular Weight Determination

Steps

Detailed Instructions

Comments



- 3 Open the default Intact Protein method and open the Deconvolute (Protein) Method Editor section.

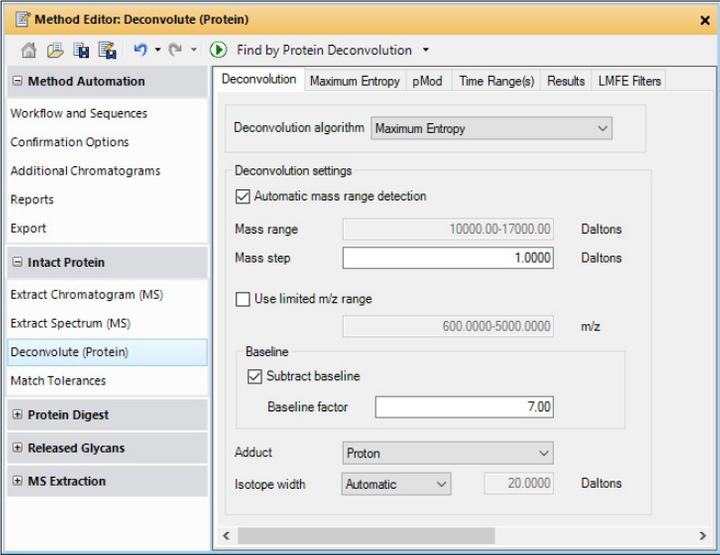
 - a Click **Method > Open**.
 - b Select *BioConfirmIntactProtein-Default.m*.
 - c Click **View > Method Editor**.
 - d Select **Intact Protein > Deconvolute (Protein)**.
 - The commands in the **View** menu toggle whether or not a window is visible. If the command is shown in blue and the button has an orange box around it, then the window is currently visible.
- 4 Select Maximum Entropy as the deconvolution algorithm.

 - On the Deconvolution tab of the Deconvolute (Protein) section of the Method Editor, verify that **Maximum Entropy** is selected for **Deconvolution algorithm**.
- 5 Verify that the Mass range is automatically detected.

 - Verify that the **Automatic mass range detection** check box is marked.
 - If you clear this check box, then you need to manually enter the Mass range which can vary for different intact proteins.

Deconvolution

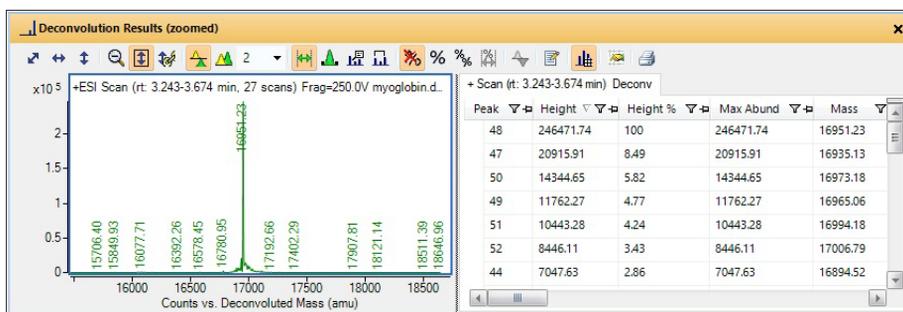
Exercise 10. Interactive Protein Molecular Weight Determination

Steps	Detailed Instructions	Comments
6 Set the Mass step to 1.	<ul style="list-style-type: none">Enter 1 for the Mass step.	
		
7 Select the extracted MS peak spectrum.	<ul style="list-style-type: none">Click the spectrum in the Biomolecule MS Spectrum window.	
8 Deconvolute the spectrum.	<ul style="list-style-type: none">Right-click the Biomolecule MS Spectrum window and click Deconvolute to start the deconvolution process.	<ul style="list-style-type: none">You can also click the arrow next to the run button in the Method Editor toolbar and select Deconvolute (Protein).
9 Review deconvolution results.	<ul style="list-style-type: none">The results appear in the Deconvolution Results window and the Biomolecules window.For information on changing the display of data in the Deconvolution Results window, see online Help.	<ul style="list-style-type: none">To compare two deconvoluted spectra, select the spectra of interest; then, click the Create Mirror Plot button,  , on the Deconvolution toolbar. If necessary, click View > Deconvolution Mirror Plot. The spectra are displayed in the Deconvolution Mirror Plot Results window. See "Exercise 11. Using the Mirror Plot window" on page 49 for more information.

Deconvolution

Exercise 10. Interactive Protein Molecular Weight Determination

Steps	Detailed Instructions	Comments
10 View peak information.	<ol style="list-style-type: none">Click the spectrum in the Deconvolution Results window to select it.Click the Spectrum Peak List button ():Click the Max Abund column heading to sort results by abundance.Click the Spectrum Peak List button () on the Deconvolution Results toolbar to close the peak list tab.	<ul style="list-style-type: none">Mass (m/z), Abundance, and Fit score are listed for each peak in the spectrum.You can change the size of the graphics pane and the table pane in the Deconvolution Results window. Select the line between them and drag it to the right or left.



11 Save the method to iii_Deconvolution_MaxEnt.m where iii are your initials	<ol style="list-style-type: none">Click Method > Save As.Enter iii_Deconvolution_MaxEnt.m for the method name.Click Save.
--	--

Exercise 11. Using the Mirror Plot window

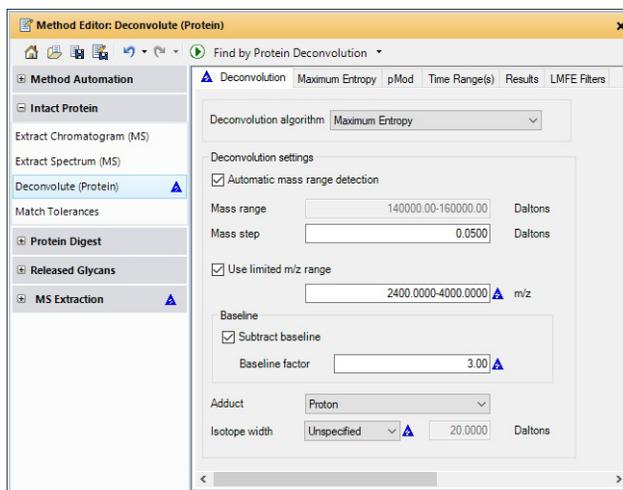
This section shows how to display a Mirror Plot of two deconvoluted biomolecules.

Steps	Detailed Instructions	Comments
1 Open the NIST mAb1.d data file.	<ul style="list-style-type: none"> a Click File > Open Data File. b Locate the NIST mAb1.d folder. c Clear Load Result Data. d Click Open. 	<ul style="list-style-type: none"> • The TIC is automatically displayed in the Sample Chromatogram Results window.
2 Open the method to use as a starting point for the new method.	<ul style="list-style-type: none"> a Click Method > Open. b Select BioConfirmIntactProtein-Default.m c Click Open. 	
3 Open the Deconvolute (Protein) Method Editor section.	<ul style="list-style-type: none"> • Select Deconvolute (Protein) from the Intact Protein section of the Method Editor. 	If the Method Editor window is not visible, click View > Method Editor to display it.

Deconvolution

Exercise 11. Using the Mirror Plot window

Steps	Detailed Instructions	Comments
<p>4 Select the deconvolution algorithm.</p> <ul style="list-style-type: none">• Use Maximum Entropy.• Use the automated mass range detection.• Use the limited m/z range of 2400 - 4000.• Use 3 for the baseline factor.	<p>a On the Deconvolution tab of the Deconvolute (Protein) section of the Method Editor, select the Deconvolution algorithm.</p> <p>b Mark the Automatic mass range detection check box.</p> <p>c Mark the Use limited m/z range check box.</p> <p>d Enter 2400 - 4000 for the <i>m/z</i> range.</p> <p>e Enter 3 for the Baseline factor.</p>	<ul style="list-style-type: none">• For more information on these parameters, press F1.
<p>5 Use the default settings for Maximum Entropy deconvolution.</p>	<ul style="list-style-type: none">• Click the Maximum Entropy tab to review settings.	
<p>6 Run the Find by Protein Deconvolution algorithm.</p>	<ul style="list-style-type: none">• Click Find and Identify > Find by Protein Deconvolution.	<p>You can also click the arrow next to the run button in the Method Editor window, and select Deconvolute (Protein).</p>



Deconvolution

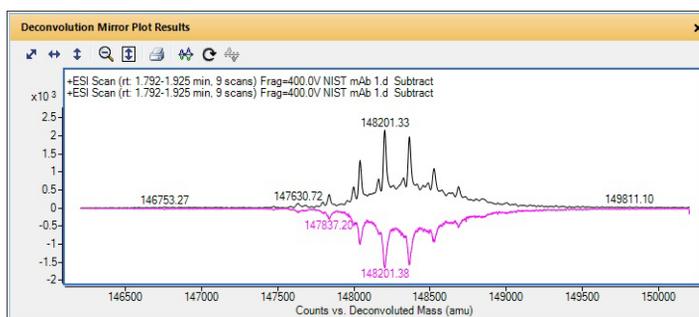
Exercise 11. Using the Mirror Plot window

Steps	Detailed Instructions	Comments
7 Review deconvolution results.	<ul style="list-style-type: none">The results appear in the Deconvolution Results window.	
8 Open the NIST mAb 2.d data file.	<ol style="list-style-type: none">Click File > Open Data File.Locate the NIST mAb 2.d sample file.Click Open.	<ul style="list-style-type: none">The TIC is automatically displayed in the Sample Chromatogram Results window.
9 Run the Find by Protein Deconvolution algorithm on NIST mAb 2.d.	<ul style="list-style-type: none">Click Find and Identify > Find by Protein Deconvolution.	
10 Review deconvolution results.	<ul style="list-style-type: none">The results appear in the Deconvolution Results window.	
11 Select both data files in the Sample Table window.	<ol style="list-style-type: none">Select one of the sample files in the Sample Table window.Press the Ctrl button and click the other sample file.	<ul style="list-style-type: none">The results for the sample files selected in the Sample Table are shown in the Deconvolution window and other windows.

Deconvolution

Exercise 11. Using the Mirror Plot window

Steps	Detailed Instructions	Comments
12 Use Mirror Plot to compare two deconvoluted spectra.	<ol style="list-style-type: none">Click the  button to show the spectra in list mode.Select a spectra from the Deconvolution window.Press the Ctrl button and select another spectra from the other data file.Click the  button to display the spectra in the Deconvolution Mirror Plot Results window.	



Exercise 12. Viewing Biomolecule Information

This exercise shows you how to view biomolecule information for deconvoluted spectra.

Steps	Detailed Instructions	Comments
1 Deconvolute myoglobin.d spectrum.	<ul style="list-style-type: none"> See “Exercise 10. Interactive Protein Molecular Weight Determination” on page 45. 	You do not need to repeat the deconvolution steps, if you have already done them in Exercise 1.
2 View the biomolecule list.	<ul style="list-style-type: none"> Click View > Biomolecules 	See Figure 5 on page 54.
3 Select the biomolecule with mass around 16973.1.	<ul style="list-style-type: none"> Click the row which has a mass around 16973.1 in the Biomolecules window. 	<ul style="list-style-type: none"> The Biomolecule MS Spectrum window and the Deconvolution Results window are both updated. A biomolecule spectrum that displays all the charge states from the original <i>m/z</i> data for that specific protein mass is shown in the Biomolecule MS Spectrum Results window.
4 Select the Biomolecule 3 spectrum in the Biomolecule MS Spectrum Results window.	<ul style="list-style-type: none"> Click the graphics area for the spectrum for Biomolecule 3. 	<ul style="list-style-type: none"> You can right-click the title of the window and click Floating. Then, you can make the window wider.
5 View the charge states found for the protein.	<ol style="list-style-type: none"> Click  on the Biomolecule MS Spectrum toolbar to show the peak information. Right-click the table and click Add/Remove Columns. Select the columns in the Available Columns list which you want to see. Click either Add or Add All ->> 	<ul style="list-style-type: none"> The following information is displayed for the ion set spectrum: <ul style="list-style-type: none"> <i>m/z</i> Abundance Charge state See Figure 6 on page 54. If you cannot see the graphics when the table is displayed, move the cursor to between the graphics and the table until it looks like . Then, click and drag to the right to increase the size of the graphics.
6 Switch from List mode to Overlay mode in the Biomolecule MS Spectrum Results window.	<ul style="list-style-type: none"> Click  on the toolbar in the Biomolecule MS Spectrum Results window. 	<ul style="list-style-type: none"> See Figure 7 on page 55.
7 Select biomolecule 1 in the biomolecule list.	<ul style="list-style-type: none"> Click the first line of the Biomolecules table. 	<ul style="list-style-type: none"> Notice that the spectrum in the Biomolecule MS Spectrum window is updated.
8 Select biomolecule 2 in the Biomolecules window.	<ul style="list-style-type: none"> Click the second line of the Biomolecules table. 	<ul style="list-style-type: none"> Notice that a different spectrum is shown in the Biomolecule MS Spectrum window.

Deconvolution

Exercise 12. Viewing Biomolecule Information

Steps	Detailed Instructions	Comments
9 Print a biomolecule report.	<ol style="list-style-type: none">Display the Reports section in the Method Editor by selecting Method Automation > Reports.Review the parameters in both the Templates and Layout tabs.Click Biomolecule Report from the File > Print menu to print the report.	<ul style="list-style-type: none">You can use either PDF-based reporting or Microsoft Excel reporting.When you print a Biomolecule Report, it uses the Intact Protein, the Protein Digest, or the Released Glycans template, depending on the workflow selected in the Sample Table window.If the workflow is Custom, then if you use the Find Peptides command, the Peptide Digest report template is used; otherwise, the Intact Protein report template is used.

Label	Mass	RT	Height	Min Z	Max Z	File	Mining Algorithm	Z Count	Use for %Quant	Fit Score
Biomolecule 1	16951.7006	3.421	4138408	6	34	myoglobin.d	Maximum Entropy D	26	<input type="checkbox"/>	10
Biomolecule 2	16934.9138	3.409	357004	6	30	myoglobin.d	Maximum Entropy D	23	<input type="checkbox"/>	9
Biomolecule 3	16974.3351	3.409	238299	6	23	myoglobin.d	Maximum Entropy D	15	<input type="checkbox"/>	9

Figure 5. Biomolecules window for myoglobin.d

Peak	m/z	Abund	Z	Max Abund	Area
1	322.048	14738.55		14707.22	777
2	652.9793	10673.62		10673.26	5261
3	663.4531	14289.77	1	14094.7	2225
4	664.4554	7362.35	1	7293.19	1808
5	678.3782	11343.33		11335.07	9083
6	679.0638	33580.73		33573.96	15789
7	679.6011	6808.4		6768.74	1494
8	685.4358	9048.16		9048.11	3312
9	693.6207	9546.3	2	9545.37	1886
10	693.8714	14461.77	2	14176.11	1823
11	694.1217	12474.57	2	12160.7	1707
12	694.3714	8719.89	2	8708.22	1388
13	700.1014	6694.05		6693.59	8414
14	704.0923	6478.82		6458.52	10830

Figure 6. Peak information for myoglobin.d displayed in the Biomolecule MS Spectrum window

Deconvolution

Exercise 12. Viewing Biomolecule Information

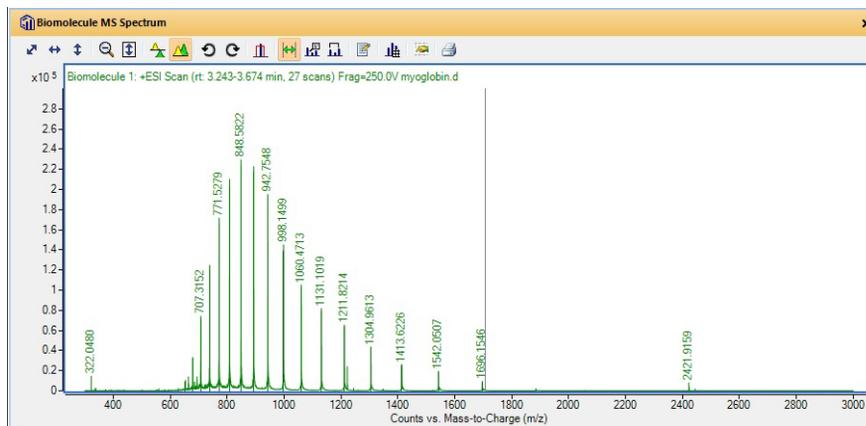


Figure 7. Biomolecule MS Spectrum Results window for myoglobin.d (Overlay Mode)

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

www.agilent.com

© Agilent Technologies, Inc. 2018

Revision A, August 2018



G6829-90028 Revision A

