

Agilent MassHunter Workstation— Unknowns Analysis

Familiarization Guide

In this Guide... 3

Before you begin these exercises 4

Task 1: Identify Compounds with TIC Analysis 5

Task 2: Identify Compounds with Deconvolution 16

Task 3: Generate the Report 39



Identify Compounds
with TIC Analysis

Identify Compounds
with Deconvolution

Generate the Report



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2015

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.

Manual Part Number

G3335-90187

Edition

January 2015

Printed in USA

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

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

If software is for use in the performance of a U.S. Government prime contract or sub-contract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14

(June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

Safety Notices

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 step-by-step exercises to help you learn to use the Unknowns Analysis program. You can do these exercises with the demonstration analysis, method, and library files, shipped with the system installation disk, or with data you acquire.

Before you begin these exercises

Copy files from the installation disk to your hard disk

1. Insert the MassHunter Quantitative Analysis installation DVD into your computer.
2. Navigate to your DVD drive:
\Data,
3. If the folder is in a compressed format, extract the data files from their zip format.
4. Copy the **Data** folder from your installation disk in uncompressed format to any location on your hard disk.

This folder contains all of the data, method, and library files needed for these exercises. Do not reuse the example data files on your system unless you know that they are identical to the originals on the disk. If the example data files already on the system do not match the original ones of the disk exactly, then the results obtained during these exercises will not match those shown in this guide.

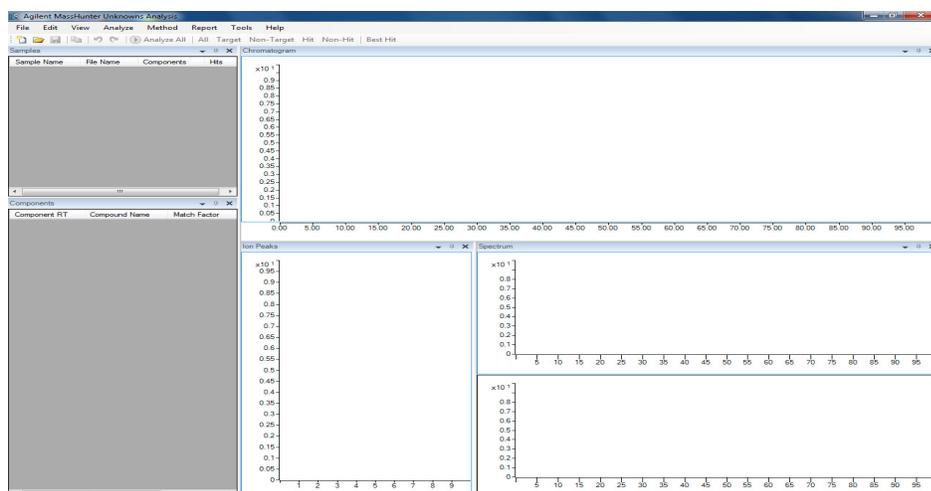
Task 1: Identify Compounds with TIC Analysis

Create a new analysis

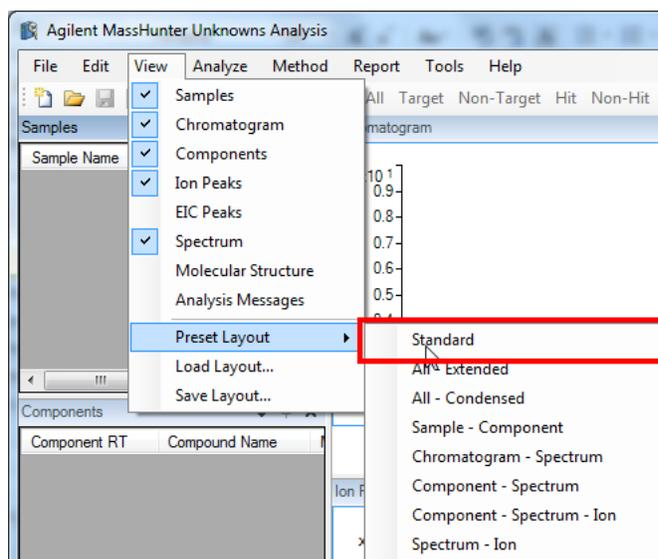
1. Start **Unknowns Analysis** by double-clicking the desktop icon.
or
Click **Start > Agilent > Quant Tools > Unknowns Analysis**.



When you open the program, the default layout appears.

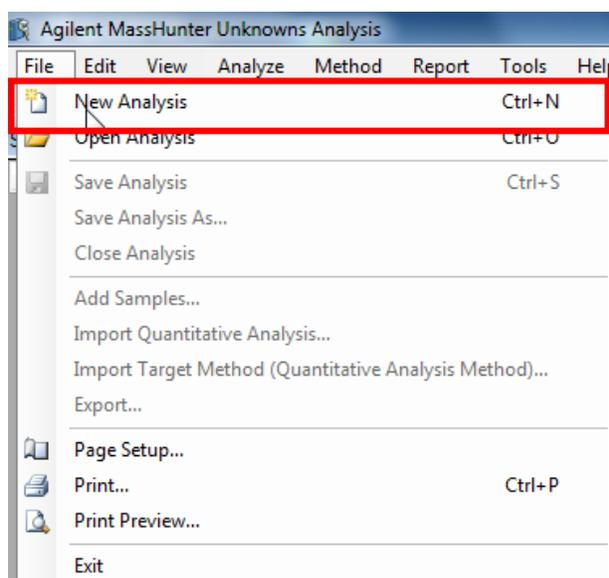


If the default layout is not present, click **View > Preset Layout > Standard** to restore the default layout before creating a new analysis.



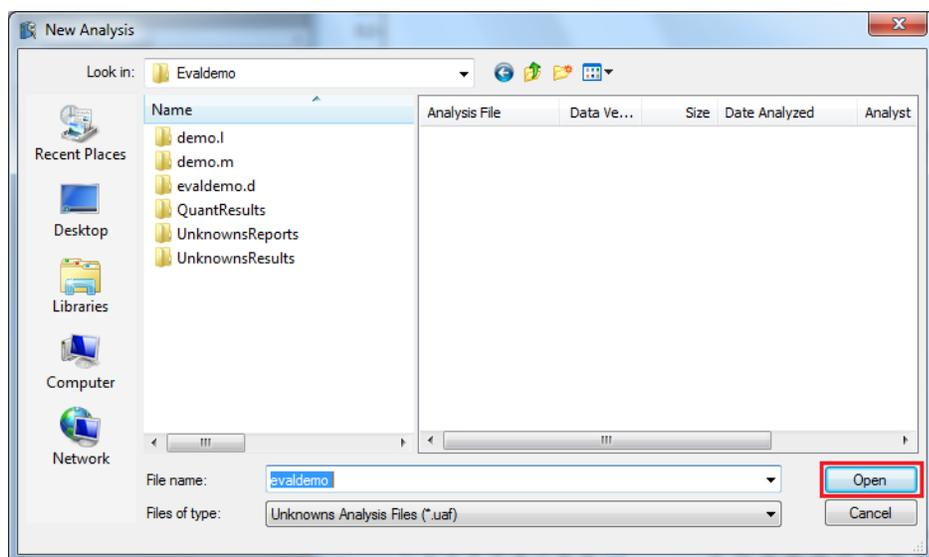
Task 1: Identify Compounds with TIC Analysis

2. Select **File > New Analysis**.



3. Navigate to **MassHunter\Data\Evaldemo**, or the folder where the data file to be analyzed is stored.

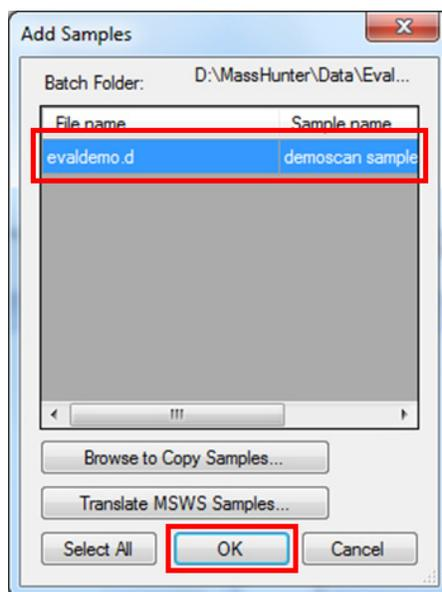
4. Type the analysis name **evaldemo** for the analysis, and click **Open**.



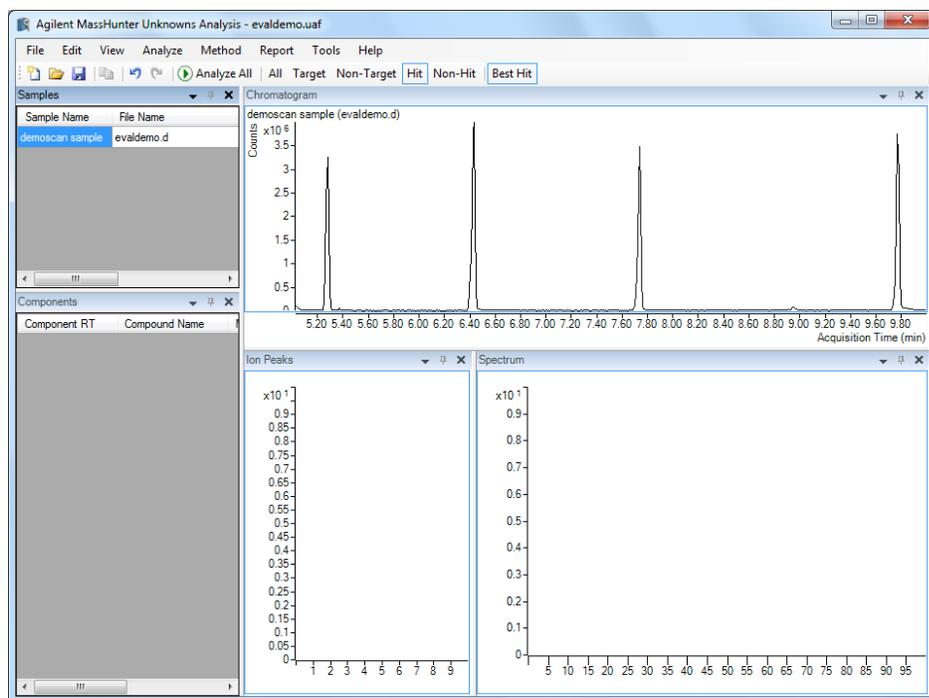
Task 1: Identify Compounds with TIC Analysis

Add samples to the analysis

1. Select **File > Add Samples**.
2. Select the sample file(s) and click **OK** to add the sample to the batch.



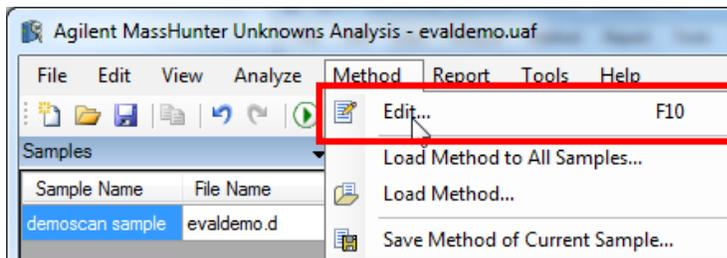
The Analysis table is no longer empty. It now contains the demo sample.



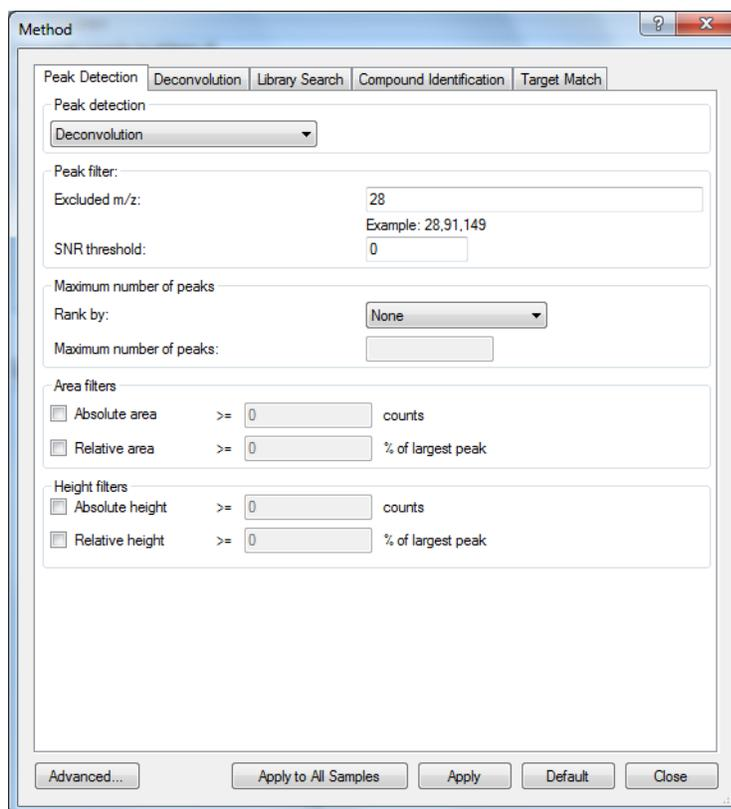
Task 1: Identify Compounds with TIC Analysis

Set up the method for the analysis

Select **Method > Edit**.



The Method dialog box standard view appears. For this task, we will use the **Standard** view.

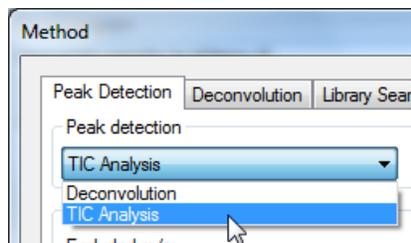


Note that these are the default parameters for the method. You can click **Default** at the bottom of the Method dialog box to restore default parameters before creating a new method in the next step.

Task 1: Identify Compounds with TIC Analysis

Set Peak Detection options

1. Select **TIC Analysis** from the **Peak detection** drop-down menu.
2. In the **Maximum number of peaks** section, select **Area** from the **Rank by** drop-down menu, and enter **5** for the **Maximum number of peaks**.
3. In the **Area filters** section, select **Relative area** and enter **1** for the **% of largest peak**.



- **TIC Analysis:** Identifies the chromatographic peaks using integration instead of deconvolution.
- **Deconvolution:** Deconvolutes the components in the chromatogram and extracts the 'clean' spectra from background noise based on both retention time and peak shape.

Maximum number of peaks

Rank by:

Maximum number of peaks:

Area filters

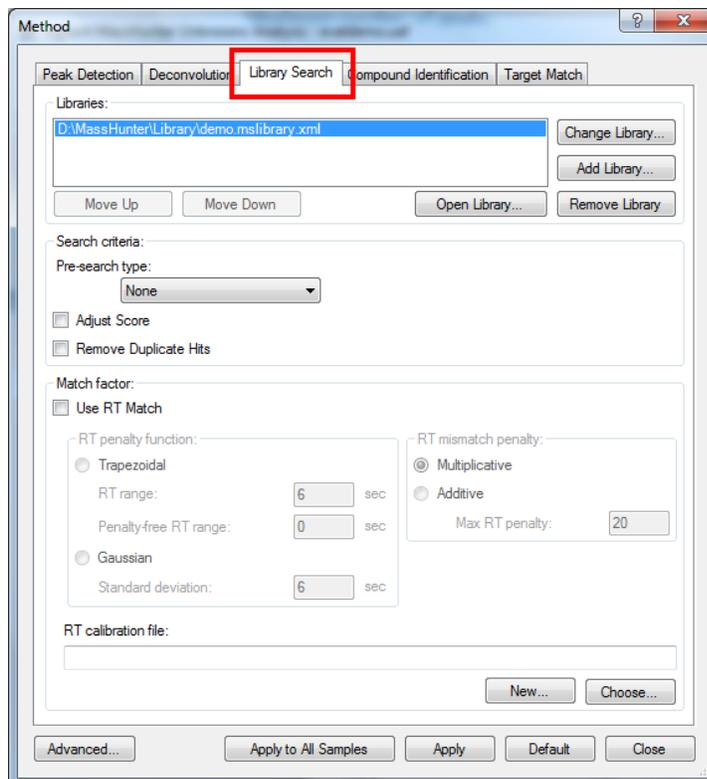
Absolute area >= counts

Relative area >= % of largest peak

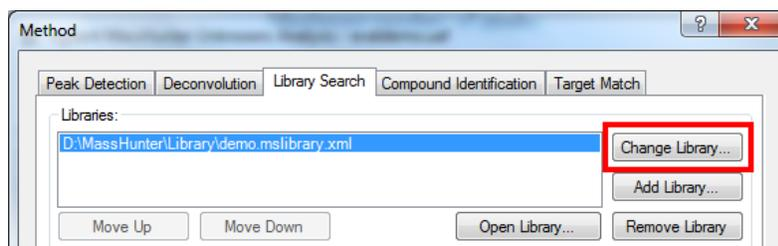
Task 1: Identify Compounds with TIC Analysis

Set Library Search options

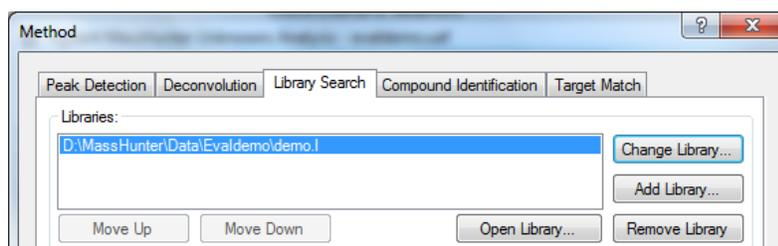
1. Click **Library Search**.



2. Click **Change Library**.



3. Navigate to **MassHunter\Data\Evaldemo**, or the relevant folder, select **demo.L**, and click **Open**.



Task 1: Identify Compounds with TIC Analysis

4. In the **Search criteria** section, select **None** from the **Pre-search** drop-down menu.

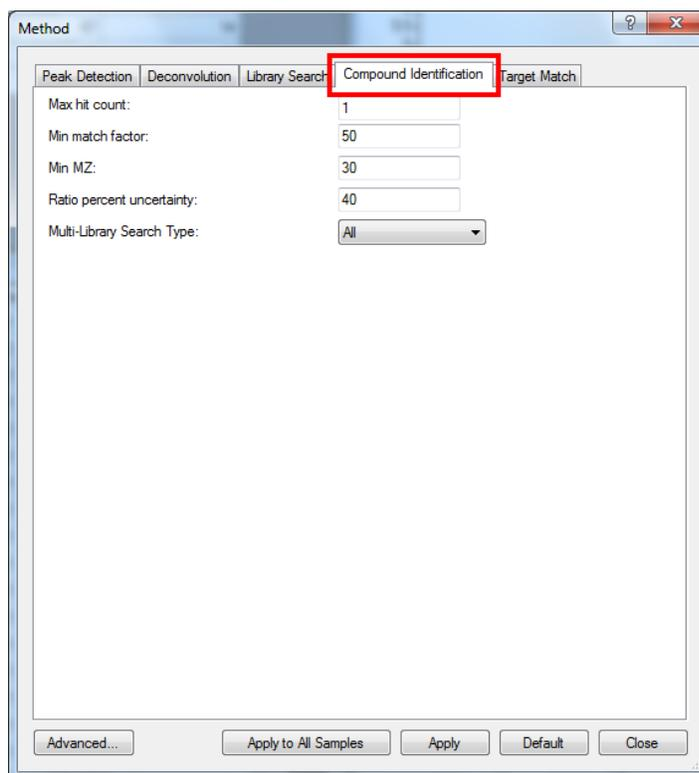


There are 3 Pre-search types: **None**, **Normal**, and **Fast**. By default, Unknowns Analysis uses **Normal**.

- **None:** The library search is not subjected to a preliminary screening process.
- **Normal:** The screening algorithm uses the entire library as the list of candidates if the indexing scheme does not produce enough candidates. It is 50-100 times faster than no pre-search, with essentially zero false negatives rate for high-scoring hits (match score > 80).
- **Fast:** The screening algorithm uses whatever list of candidates it gets from the index and avoids the entire library-search even if there are not enough candidates found. The speed is 100-1000 times faster than no pre-search, with $\geq 1\%$ false negatives rate for high-scoring hits.

Set Compound Identification options

1. Click **Compound Identification**.



For this task, we will use the default Compound Identification parameters.

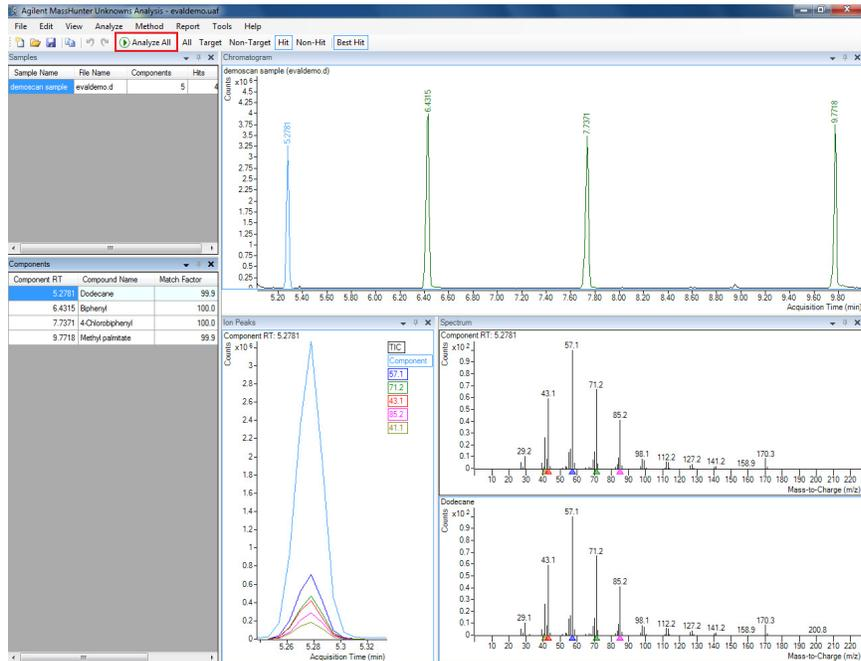
2. Click **Apply to All Samples**, and then click **Close**.

Task 1: Identify Compounds with TIC Analysis

Analyze and review results

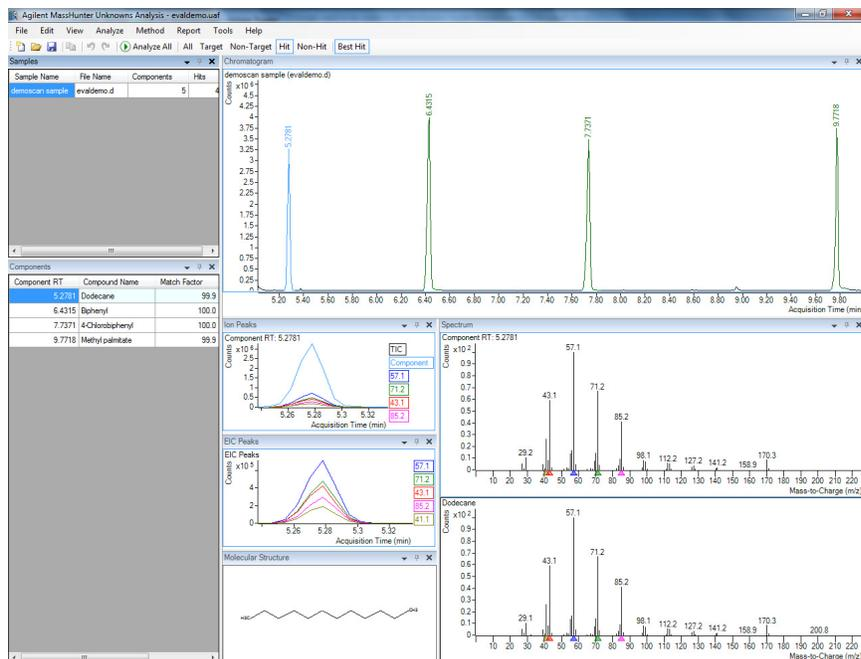
1. Click **Analyze All**.

After the analysis is complete, the main view that appears should look like the example below. This is the default layout and contains the default column settings. If you see a different layout than the one in the example below, select **View > Preset Layout > Standard** to reset the standard layout.



2. Select **View > Preset Layout > All-condensed**.

The system displays the All-condensed view.



Task 1: Identify Compounds with TIC Analysis

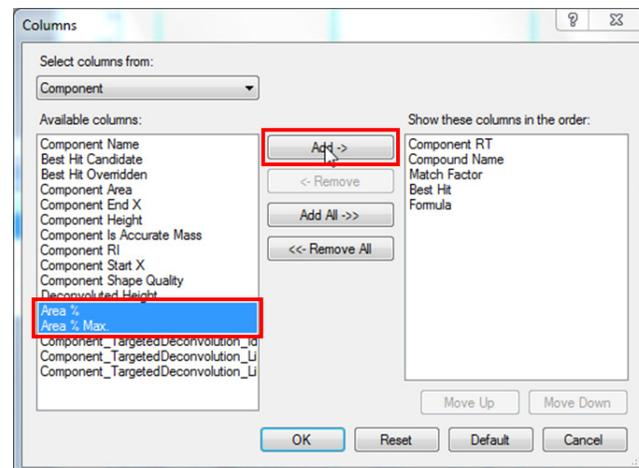
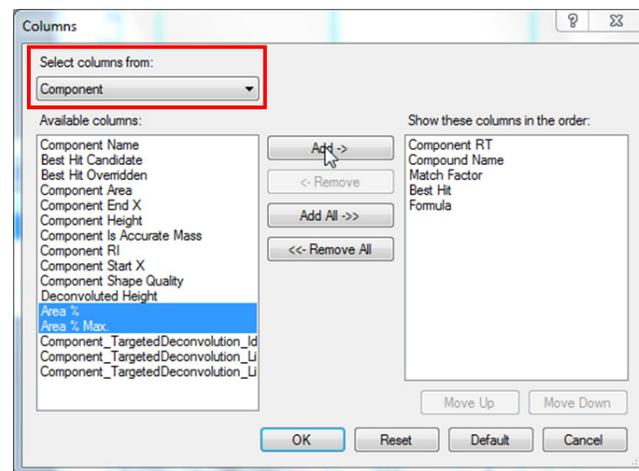
3. Select the **demSCAN sample** from the **Sample** table.
4. Right-click any column header in the **Components** window, and select **Add/Remove Columns**.
5. Select **Component** from the **Select columns from** drop-down menu.
6. Select **Area %** and **Area % Max** from the **Available columns** list, and click **Add**.

Click one of the following toolbar buttons to view the changes in the **Components** window:

- **All**: View all the peaks.
- **Hit**: View the peaks that are found in the library search.
- **Non-Hit**: View the peaks that are not found in the library search.

The screenshot shows the 'Components' window with a table of peaks. A right-click context menu is open over the 'Match Factor' column header, with 'Add/Remove Columns...' highlighted.

Component RT	Compound Name	Match Factor
5.2781	Dodecane	9
6.4315	Biphenyl	10
7.7371	4-Chlorobiphenyl	10
9.7718	Methyl palmitate	9



- **Area %**: Percentage of the peak area sum
- **Area % Max**: Percentage of the largest peak area

Task 1: Identify Compounds with TIC Analysis

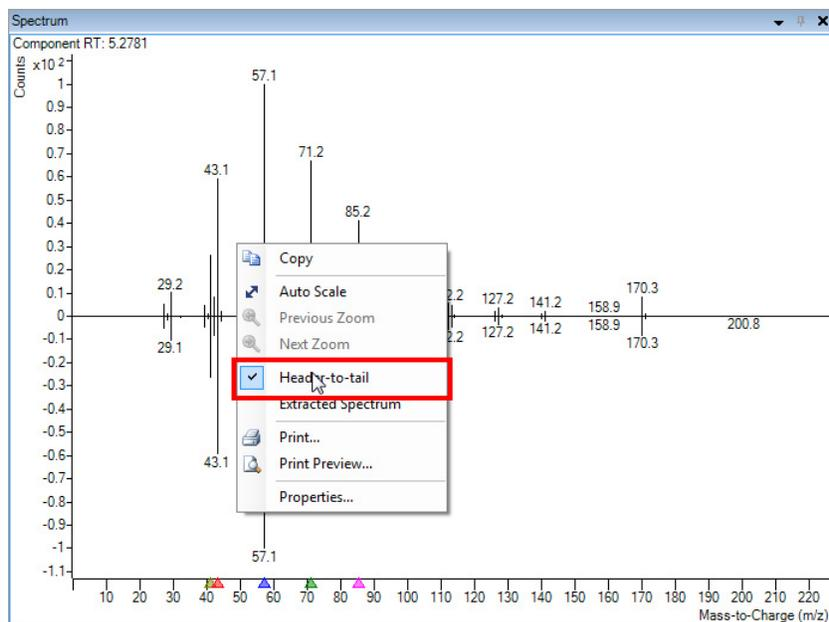
- Verify that the selected columns are moved to the **Show these columns in the order** list, and click **OK**.
- From the **Components** table, select a component in the **Component RT** column.

Component RT	Compound Name	Match Factor	Best Hit	Formula	Area %	Area % Max.
5.2781	Dodecane	99.9	<input checked="" type="checkbox"/>	C ₁₂ H ₂₆	19.158	63.44
6.4315	Biphenyl	100.0	<input checked="" type="checkbox"/>	C ₁₂ H ₁₀	30.196	100.00
7.7371	4-Chlorobiphenyl	100.0	<input checked="" type="checkbox"/>	C ₁₂ H ₉ Cl	25.572	84.69
9.7718	Methyl palmitate	99.9	<input checked="" type="checkbox"/>	C ₁₇ H ₃₄ O ₂	23.731	78.59

View the **Chromatogram**, **Spectrum**, **Ion Peaks**, **EIC Peaks**, and **Molecular Structure** for the selected component.

In the **Spectrum** window, the top spectrum is from the component, and the bottom spectrum is from the library. The **Match Factor** in the **Components** table reflects how closely the two spectrum match.

To change to the Header-to-tail view, right-click inside the **Spectrum** window and select **Header-to-tail**.



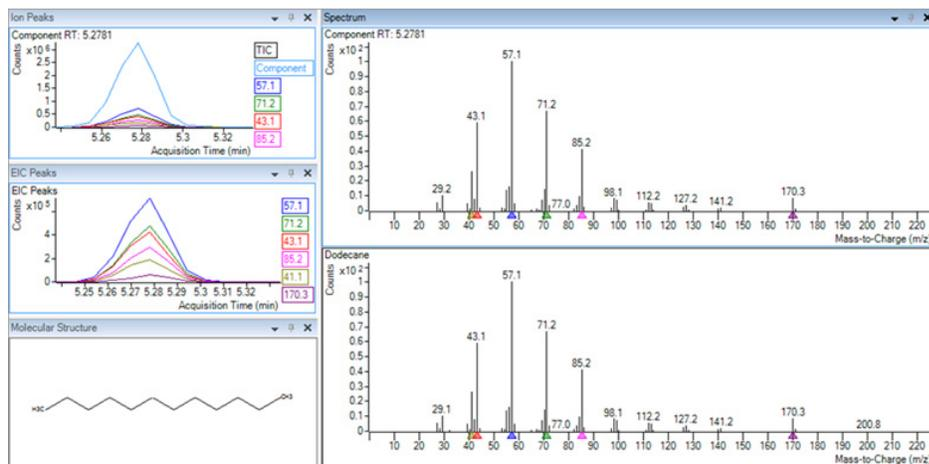
The **Ion Peaks** and **EIC Peaks** windows show the extracted chromatograms of the selected ions. The EIC traces and their numeric identifiers to the right of the display are color-coded.

To interactively add the ion chromatogram traces in the **Ion Peaks** and **EIC Peaks** window to the display, click in any **Mass Spectral Display** area of the **Spectrum** window. If the selected m/z chromatogram is not already displayed, it will be added to the **Ion Peaks** and **EIC Peaks** window and a ▲ symbol of the same color will be at the appropriate m/z position below the x-axis in the **Spectrum** window.

Task 1: Identify Compounds with TIC Analysis

To remove an ion chromatogram trace (and its numeric identifier) from the **Ion Peaks** and **EIC Peaks** window, click on its numeric identifier or on the corresponding m/z value position in the **Spectrum** window.

The **Molecular Structure** is from the library. If the searched library does not contain the structures for the entries, nothing will be displayed in the **Molecular Structure** window.



9. To save the analysis, select **File > Save Analysis**.

10. Click **File > Exit**.

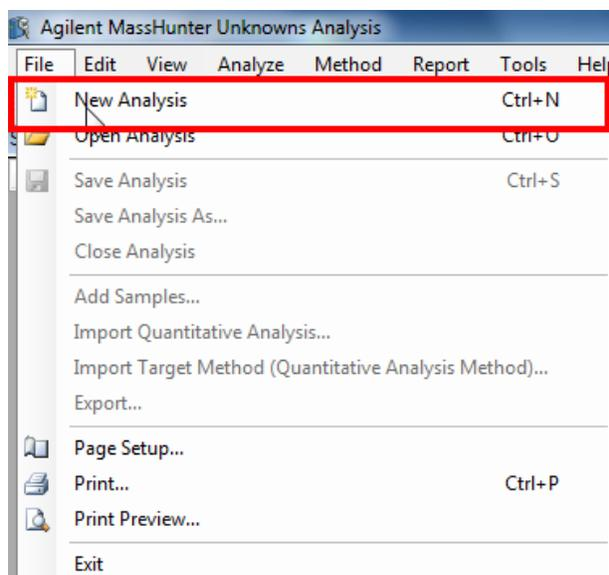
Task 2: Identify Compounds with Deconvolution

Create a new analysis

1. Start **Unknowns Analysis** by double-clicking the desktop icon.
or
Click **Start > Agilent > Quant Tools > Unknowns Analysis**.
2. Select **File > New Analysis**.



Unknowns Analysis



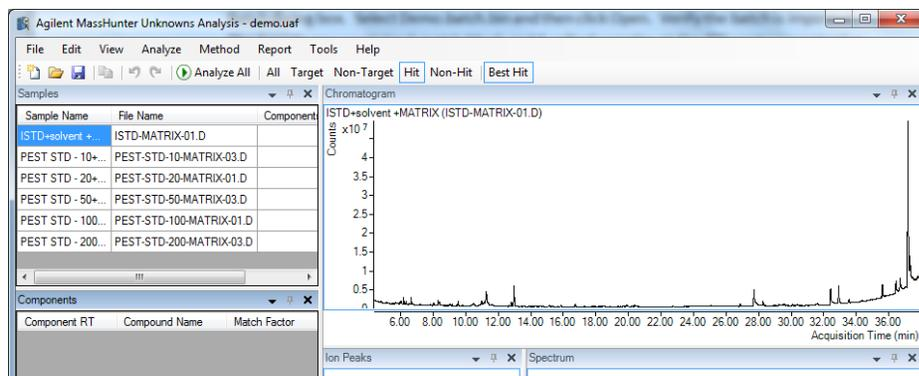
3. Navigate to **\Your Directory\
RI-PEST-MATRIX**.
4. Type the analysis name **demo**,
and click **Open**.

Task 2: Identify Compounds with Deconvolution

Add samples to the analysis

1. Select **File > Import Quantitative Analysis**.
2. Select **Demo.batch.bin**, and click **Open**.

Verify the batch is imported. The **Sample** window now contains one matrix blank and five spiked samples at the different concentration levels. The **Chromatogram** shows the TIC of the sample selected in the **Sample** window.



Set up the method for the analysis

Press **F10** or select **Method > Edit**.

Task 2: Identify Compounds with Deconvolution

Set Peak Detection options

Select **Deconvolution** from the **Peak detection** drop-down menu, and click **Apply to All Samples**.

The screenshot shows the 'Method' dialog box with the 'Peak Detection' tab selected. The 'Peak detection' dropdown menu is set to 'Deconvolution' and is highlighted with a red box. Below this, there are several input fields: 'Excluded m/z' (28), 'SNR threshold' (0), 'Maximum number of peaks' (None), 'Rank by' (None), and 'Maximum number of peaks' (empty). There are also checkboxes for 'Area filters' (Absolute area, Relative area) and 'Height filters' (Absolute height, Relative height), each with a '>=' input field set to 0. At the bottom, the 'Apply to All Samples' button is highlighted with a red box.

Set Deconvolution options

1. Click **Deconvolution**.

The screenshot shows the 'Method' dialog box with the 'Deconvolution' tab selected. The 'Deconvolution' tab is highlighted with a red box. The 'Resolution' field is empty. The 'RT window size factor' is 25,50,100,200. The 'Extraction window' section includes 'Left m/z delta' (0.3), 'Right m/z delta' (0.7), and 'm/z delta units' (AMU). There are checkboxes for 'Use integer m/z values' and 'Use base peak shape'. The 'Sharpness threshold' is 25%. The 'Ion peaks' section includes 'Min # of ion peaks' (3) and 'Max # of ion peak shapes to store' (10). At the bottom, the 'Apply to All Samples' button is highlighted with a red box.

Task 2: Identify Compounds with Deconvolution

The default parameters for deconvolution display. By default, there are four values (25, 50, 100, 200) for the **RT window size factor**. Select any set of Window Size Factor (WSF) values in a comma-separated format.

The screenshot shows the 'Method' dialog box with the 'Deconvolution' tab selected. The 'Resolution' section is highlighted with a red box, showing the 'RT window size factor' set to '25,50,100,200'. Other parameters include 'Extraction window' (Left m/z delta: 0.3, Right m/z delta: 0.7, m/z delta units: AMU), 'Component shape' (Use base peak shape: unchecked, Sharpness threshold: 25%), and 'Ion peaks' (Min # of ion peaks: 3, Max # of ion peak shapes to store: 10). Buttons at the bottom include 'Advanced...', 'Apply to All Samples', 'Apply to Selected Sample', 'Default', and 'Close'.

2. In the **Extraction Window** section, select **Use integer m/z values**.

The WSF represents a dimensionless scale of the correlation window for grouping ion peaks into components, equivalent to Resolution and AMDIS. A smaller value (higher resolution) separates closely spaced peaks, finds more components, and runs longer. A larger value is used for wider peaks. Using multiple values covers all kinds of peaks without manual optimization.

This close-up shows the 'Extraction window' section with 'Left m/z delta' at 0.3, 'Right m/z delta' at 0.7, and 'm/z delta units' set to 'AMU'. The checkbox for 'Use integer m/z values' is checked and highlighted with a red box.

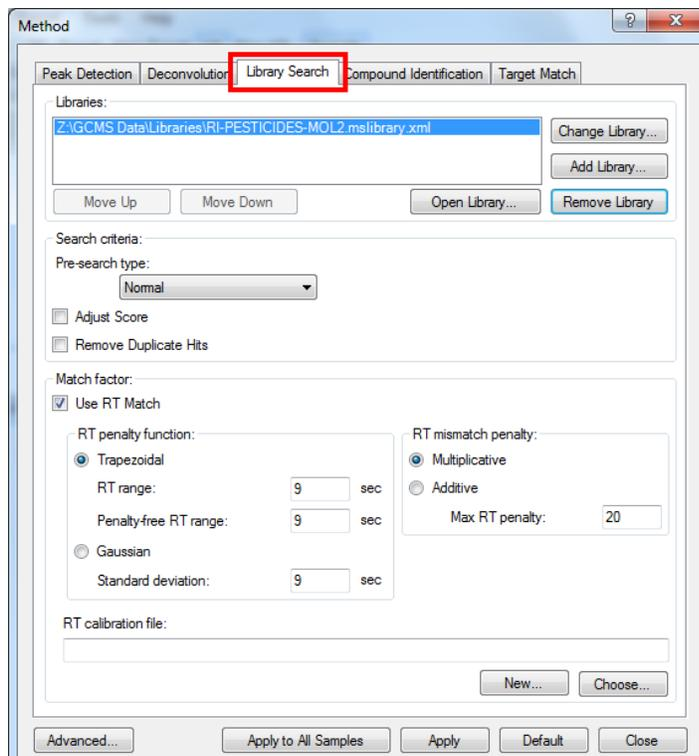
3. Click **Apply to All Samples**.

Use integer m/z values runs the deconvolution with both integer and filtered m/z , and provides the best results.

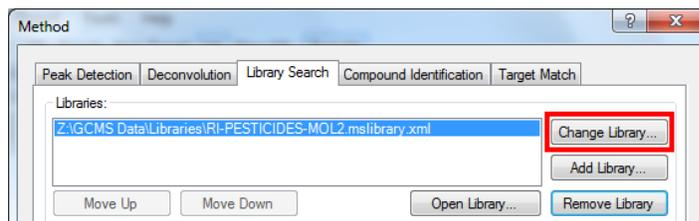
Task 2: Identify Compounds with Deconvolution

Set Library Search options

1. Click **Library Search**.



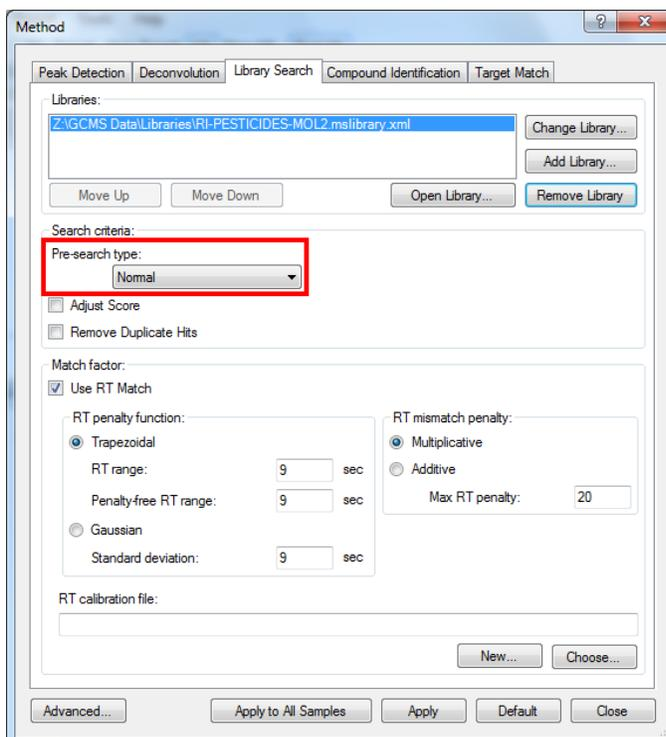
2. Click **Change Library**.



3. Navigate to the relevant folder, select **RI-PESTICIDES-MOL2.mslibrary.xml**, and click **Open**.

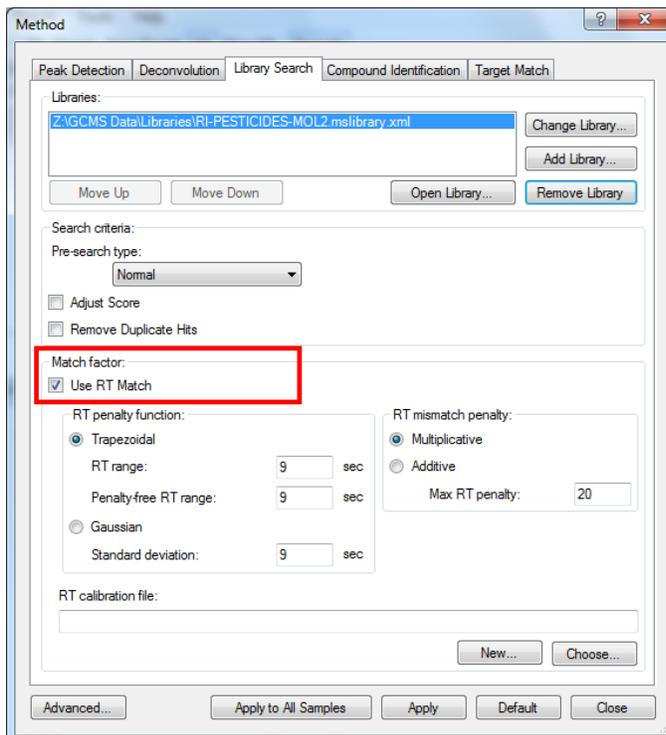
Task 2: Identify Compounds with Deconvolution

4. In the **Search criteria** section, select **Normal** from the **Pre-search type** drop-down menu.



- Select **Adjust Score** to give the closest library match scores to NIST.
- Select **Remove Duplicate Hits** to remove duplicate hits that appear in the hit list for a given target spectrum. This deals with duplicate and highly similar library entries such as seen in NIST, and only returns the single library entry with the highest fit score.

5. In the **Match factor** section, select **Use RT Match**.



Task 2: Identify Compounds with Deconvolution

- In the **RT penalty function** section, select **Trapezoidal** and enter the following:

RT range: 9

Penalty-free RT range: 9

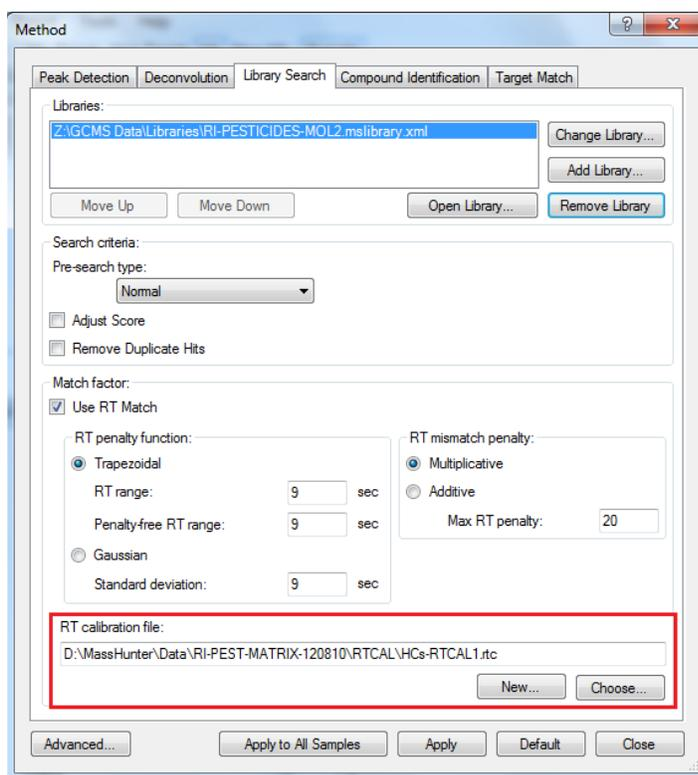
The screenshot shows the 'Method' dialog box with the 'Deconvolution' tab selected. The 'Libraries' section shows a list with 'Z:\GCMS Data\Libraries\RI-PESTICIDES-MOL2.mslibrary.xml'. The 'Search criteria' section has 'Pre-search type' set to 'Normal'. The 'Match factor' section has 'Use RT Match' checked. Under 'RT penalty function', the 'Trapezoidal' radio button is selected and highlighted with a red box. The 'RT range' is set to 9 sec and the 'Penalty-free RT range' is also set to 9 sec. The 'RT mismatch penalty' section has 'Multiplicative' selected and 'Max RT penalty' set to 20. The 'RT calibration file' section is empty. At the bottom, the 'Choose...' button is visible.

- In the RT calibration file section, click **Choose**.

This screenshot is identical to the one above, showing the 'Method' dialog box with the 'Deconvolution' tab. The 'RT penalty function' section is still highlighted with a red box. In this view, the 'Choose...' button in the 'RT calibration file' section is highlighted with a red box, indicating the next step in the process.

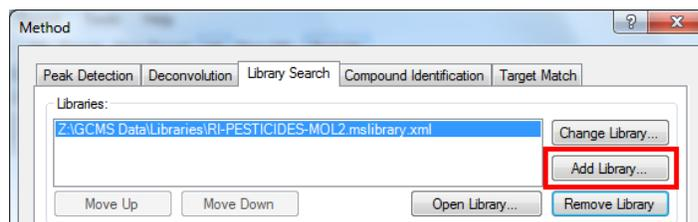
Task 2: Identify Compounds with Deconvolution

8. Navigate to the relevant folder, and select **HCS-RTCAL1.rtc**.



RT/RI calculation is used with library matching to lower the false positive rate. The window is set to ± 9 seconds to qualify the hits from the Library Search.

9. In the **Libraries** section, click **Add Library**.

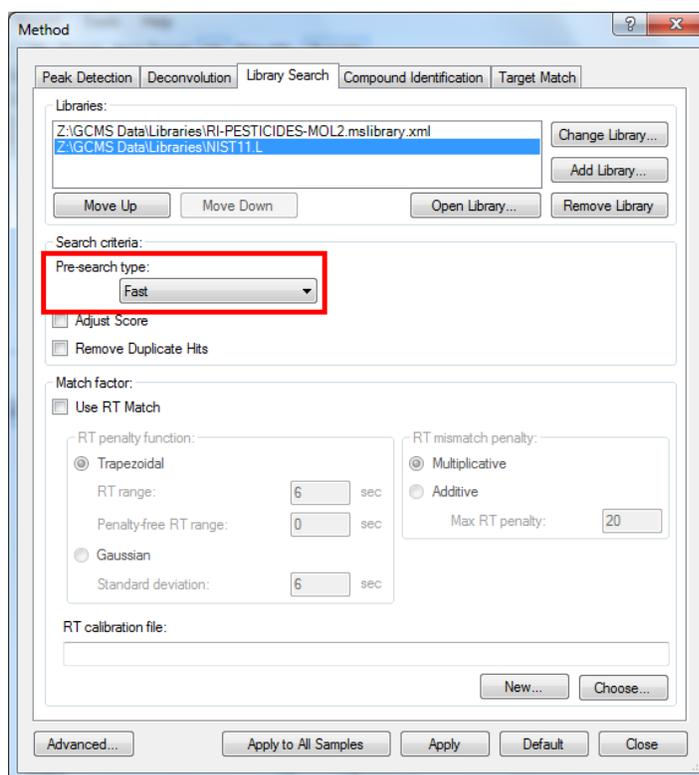


10. Navigate to the relevant folder, and select **NIST11.L**.

Multiple libraries can be used in Library Search. For this example, the target MS library contains 900+ pesticides with Retention Indexes (RI) information. NIST11.L can be used for the additional confirmation.

Task 2: Identify Compounds with Deconvolution

11. Select **Fast** from the **Pre-search type** drop-down menu.



12. Click **Apply to All Samples**.

You are able to set different Library Search parameters for different libraries.

Task 2: Identify Compounds with Deconvolution

Set Compound Identification options

Click **Compound Identification**.

The screenshot shows a software window titled 'Method' with several tabs: 'Peak Detection', 'Deconvolution', 'Library Search', 'Compound Identification', and 'Target Match'. The 'Compound Identification' tab is selected and highlighted with a red box. The settings for this tab are as follows:

Parameter	Value
Max hit count:	1
Min match factor:	50
Min MZ:	30
Ratio percent uncertainty:	40
Multi-Library Search Type:	All

At the bottom of the dialog, there are five buttons: 'Advanced...', 'Apply to All Samples', 'Apply', 'Default', and 'Close'.

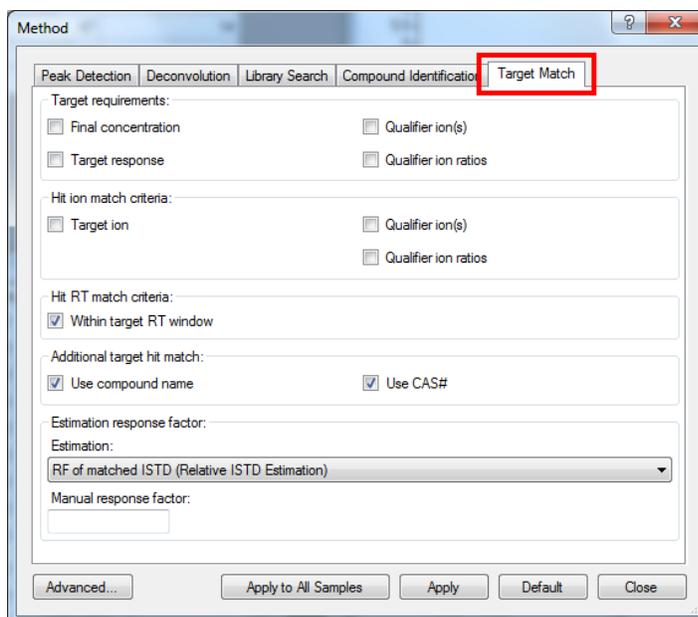
For this example, the **Min match factor** is set to 50 for the compound identification from the Library Search.

- **Max hit count:** The maximum number of Library Search hits to report per component.
- **Min MZ:** The lower m/z limit for library match score calculation.
- **Ratio percent uncertainty:** Only applicable when **Pre-search type** is selected in Library Search. The larger the value, the more Library Search candidates are generated, and the longer the library search process.
- **All:** Search all libraries (default)
- **Multi-Library Search Type:** If multiple libraries were used, two search modes are available:
 - **All:** Search all libraries (default)
 - **StopWhenFound:** Stop searching the library when enough candidates are found

Task 2: Identify Compounds with Deconvolution

Set Target Match options

1. Click **Target Match**.



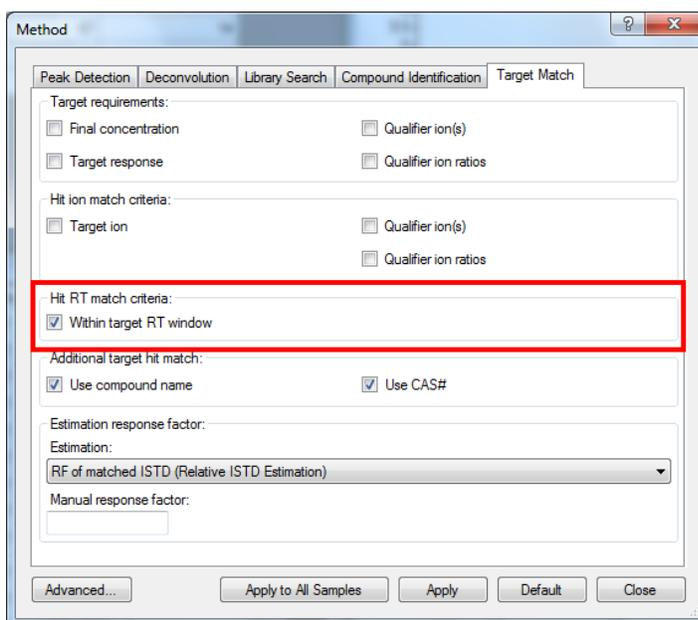
The screenshot shows the 'Method' dialog box with the 'Target Match' tab selected. The 'Target Match' tab is highlighted with a red box. The dialog contains several sections with checkboxes and a dropdown menu:

- Target requirements:**
 - Final concentration
 - Target response
 - Qualifier ion(s)
 - Qualifier ion ratios
- Hit ion match criteria:**
 - Target ion
 - Qualifier ion(s)
 - Qualifier ion ratios
- Hit RT match criteria:**
 - Within target RT window
- Additional target hit match:**
 - Use compound name
 - Use CAS#
- Estimation response factor:**
 - Estimation: RF of matched ISTD (Relative ISTD Estimation)
 - Manual response factor: []

Buttons at the bottom: Advanced..., Apply to All Samples, Apply, Default, Close.

Target Match identifies quantitation targets using the quantitation method. The goal of identifying non-target compounds is simplified by filtering out the target matches. RT window, compound name, and CAS# can be applied for **Target Match**.

2. In the **Hit RT match criteria** section, select **Within target RT window**.



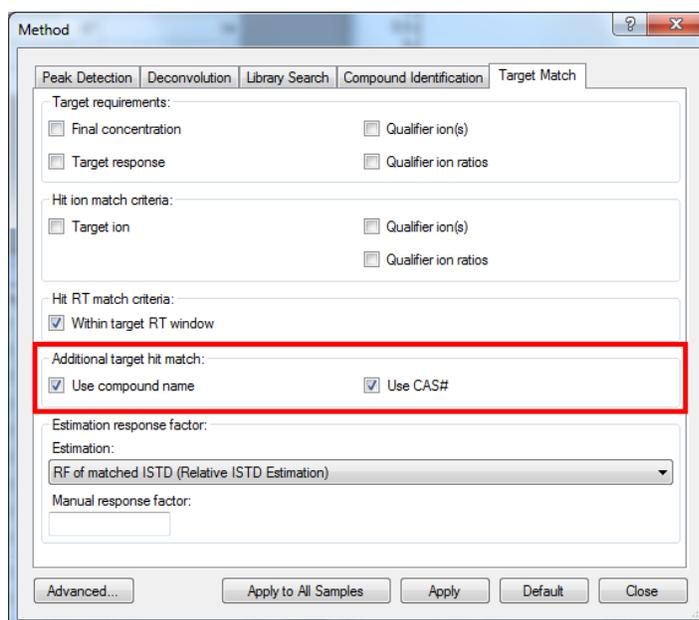
This screenshot is similar to the previous one, but the 'Within target RT window' checkbox in the 'Hit RT match criteria' section is now checked and highlighted with a red box. The 'Target Match' tab is also highlighted with a red box.

- Hit RT match criteria:**
 - Within target RT window

Buttons at the bottom: Advanced..., Apply to All Samples, Apply, Default, Close.

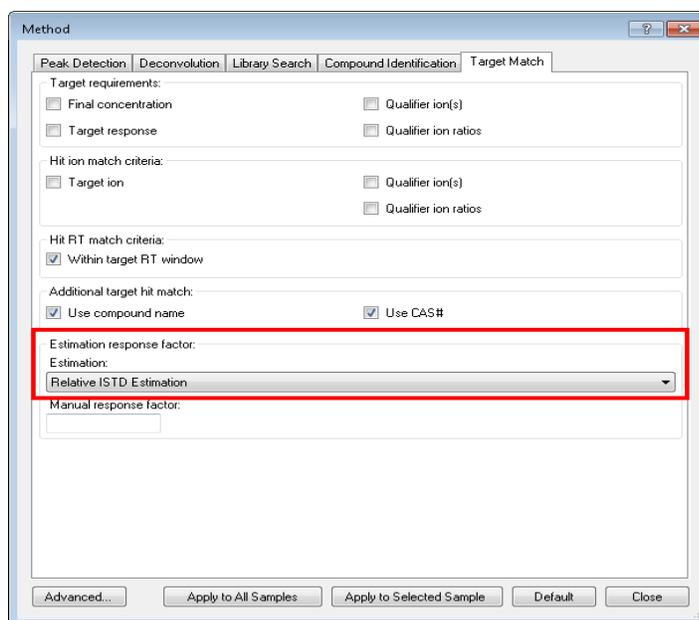
Task 2: Identify Compounds with Deconvolution

3. In the **Additional target hit match** section, select **Use compound name** and **Use CAS#**.



The screenshot shows the 'Method' dialog box with the 'Target Match' tab selected. The 'Additional target hit match' section is highlighted with a red box. In this section, the checkboxes for 'Use compound name' and 'Use CAS#' are both checked. Other sections include 'Target requirements', 'Hit ion match criteria', 'Hit RT match criteria', and 'Estimation response factor'.

4. In the **Estimation response factor** section, select **RF of matched ISTD** from the **Estimation** drop-down menu.

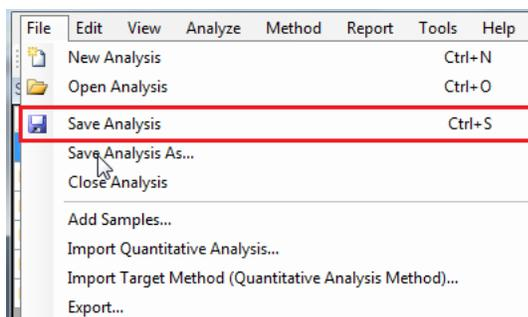


The screenshot shows the 'Method' dialog box with the 'Target Match' tab selected. The 'Estimation response factor' section is highlighted with a red box. In this section, the 'Estimation' drop-down menu is set to 'Relative ISTD Estimation'. Other sections include 'Target requirements', 'Hit ion match criteria', 'Hit RT match criteria', and 'Additional target hit match'.

Concentration estimation leverages the Quant target **Response Factors (RF)**, which are applied to Non-Target hits as well. Estimation of response factors is flexible, and can be adjusted to suit the particular analytical requirements.

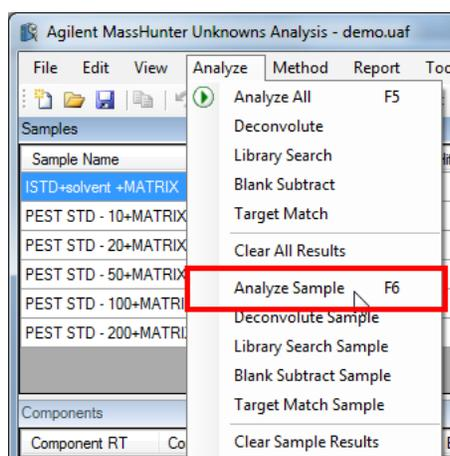
Task 2: Identify Compounds with Deconvolution

5. Click **Apply to All Samples**, and then click **Close**.
6. To save the analysis, select **File > Save Analysis**.

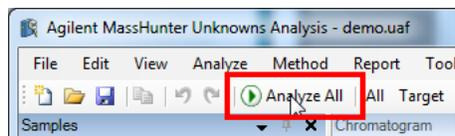


Analyze and review results

1. In the **Sample** window, select the sample **ISTD+solvent+MATRIX**.
2. Select **Analyze > Analyze Sample**.



To analyze the rest of the sample, click **Analyze All**. The analysis starts from where it left off and skips the sample(s) previously analyzed if no parameter in the method has been changed.

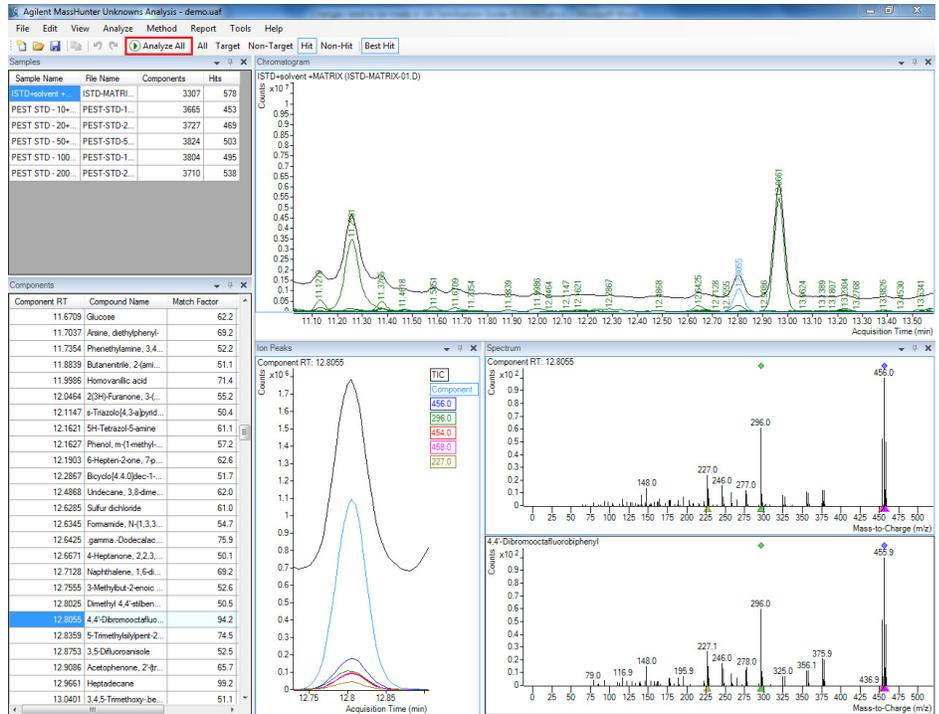


Task 2: Identify Compounds with Deconvolution

View validation information in the Analysis Messages.

Type	Target	Message
	Sample ISTD+solvent +MATRIX	Deconvolution process has been already performed before. Skipping deconvolution process.
	Sample ISTD+solvent +MATRIX	Library search process has been already performed before. Skipping library search process.
	Sample ISTD+solvent +MATRIX	Blank subtraction process has been already performed before. Skipping blank subtraction process.
	Sample ISTD+solvent +MATRIX	Target match process has been already performed before. Skipping target match process.

After the analysis is complete, the main view that appears should look like the example below. This is the default layout and contains the default column settings.



Review best hit results

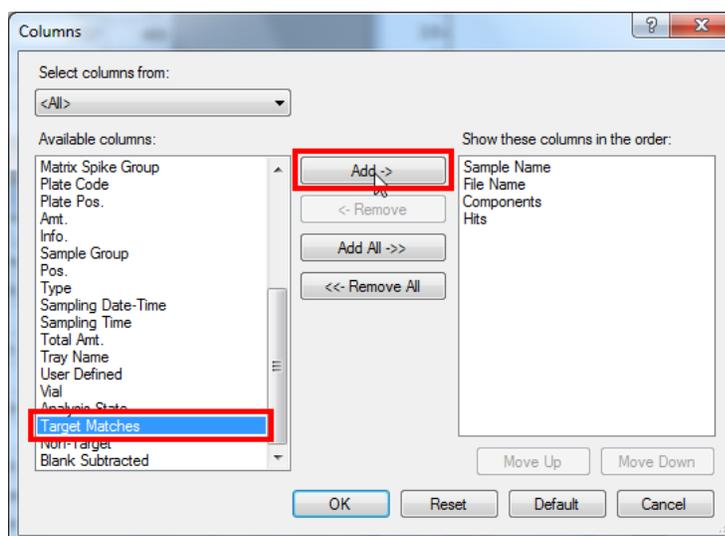
1. Right-click any column header in the **Samples** window, and select **Add/Remove Columns**.

The screenshot shows the **Samples** window with a context menu open over the **Components** column header. The menu options are:

- Add/Remove Columns...
- Analyze Sample
- Deconvolute Sample
- Library Search Sample
- Blank Subtract Sample
- Target Match Sample
- Clear Sample Results
- Delete Sample
- Format Column...
- Print...
- Print Preview...

Task 2: Identify Compounds with Deconvolution

2. Select **Target Matches** from the **Available columns** list and click **Add**.



3. Verify that the selected column is moved to the **Show these columns in the order** list, and click **OK**.

The **Target Matches** column is added to the **Samples** window.

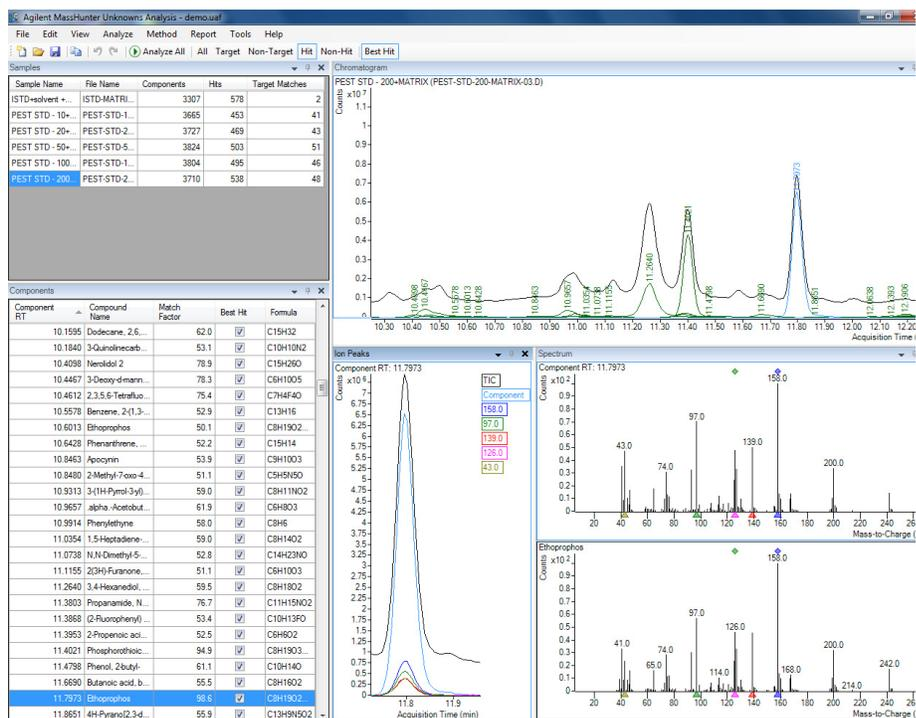
Sample Name	File Name	Components	Hits	Target Matches
ISTD+solvent +...	ISTD-MATRI...	3307	578	2
PEST STD - 10+...	PEST-STD-1...	3665	453	41
PEST STD - 20+...	PEST-STD-2...	3727	463	43
PEST STD - 50+...	PEST-STD-5...	3824	503	51
PEST STD - 100...	PEST-STD-1...	3804	493	46
PEST STD - 200...	PEST-STD-2...	3710	533	48

Task 2: Identify Compounds with Deconvolution

- Select the last sample in the **Samples** window.

Click one of the following toolbar buttons to view the changes in the **Components**, **Chromatogram**, **Ion Peaks**, and **Spectrum** windows:

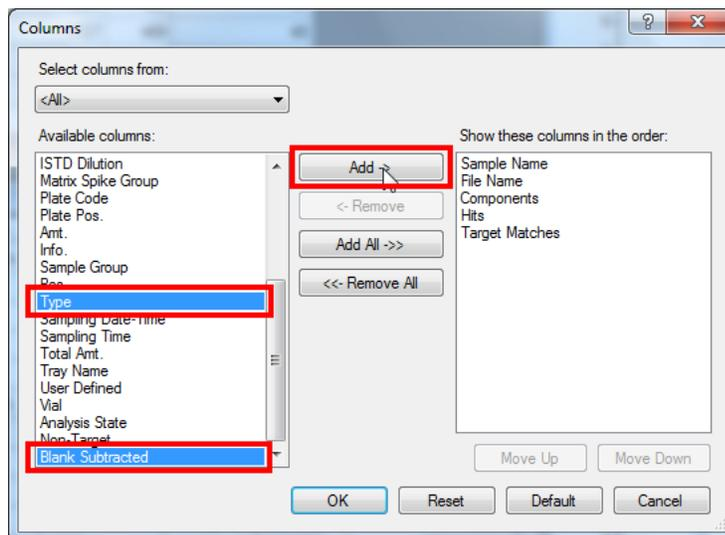
- All:** View all the peaks.
- Target:** View the peaks that are also in the quantitation method.
- Non-Target:** View the peaks that are not in the quantitation method.
- Hit:** View the peaks that are found in the library search.
- Non-Hit:** View the peaks that are not found in the library search.
- Best Hit:** View the component with the highest library match score among the multiple hits of the same compound from different resolutions.



Task 2: Identify Compounds with Deconvolution

Review blank hit subtraction results

1. Right-click any column header in the **Samples** window, and select **Add/Remove Columns**.
2. Select **Type** and **Blank Subtracted** from the **Available columns** list, and click **Add**.

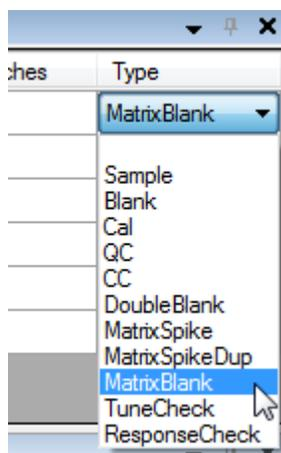


3. Verify that the selected columns are moved to the **Show these columns in the order** list, and click **OK**.

The **Type** and **Blank Subtracted** columns are added to the **Samples** window.

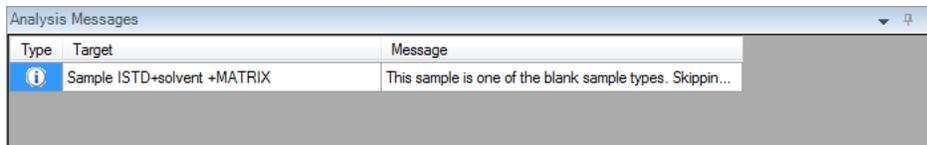
Sample Name	File Name	Components	Hits	Target Matches	Type	Blank Subtracted
ISTD+solvent +MATRIX	ISTD-MATRI...	3307	578	2	MatrixBlank	0
PEST STD - 10+MATRIX	PEST-STD-1...	3665	453	41	Cal	183
PEST STD - 20+MATRIX	PEST-STD-2...	3727	469	43	Cal	167
PEST STD - 50+MATRIX	PEST-STD-5...	3824	503	51	Cal	153
PEST STD - 100+MATRIX	PEST-STD-1...	3804	495	46	Cal	160
PEST STD - 200+MATRIX	PEST-STD-2...	3710	538	48	Cal	149

4. Note the list of available samples in the **Type** drop-down menu.



Task 2: Identify Compounds with Deconvolution

The values shown in the **Blank Subtracted** column in the **Samples** window represent the number of hits that were blank subtracted from the samples. Verify that the message “This sample is one of the blank sample types. Skipping blank subtraction process.” for the **ISTD+solvent+MATRIX** sample appears in the **Analysis Messages** window.



Type	Target	Message
i	Sample ISTD+solvent +MATRIX	This sample is one of the blank sample types. Skippin...

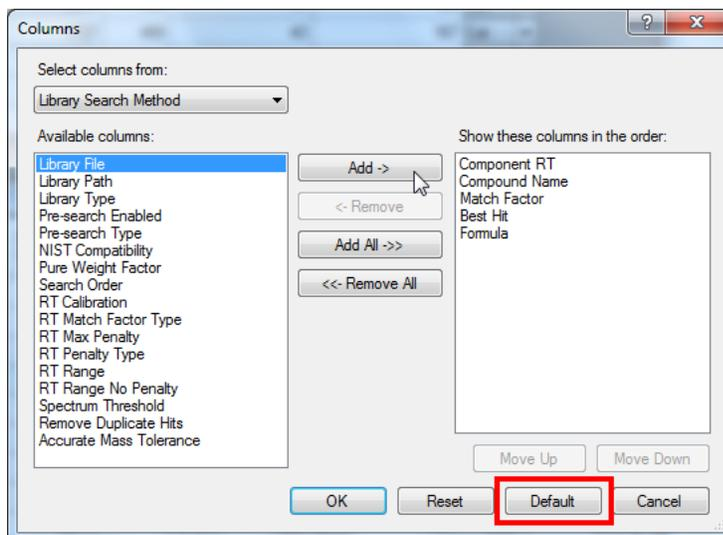
Close the **Analysis Messages** window.

Hits in a sample are marked as **Blank Subtracted Hits** when the same hit is found in the blank with $RT \pm 5FWHM$. FWHM of a typical GC-MS peak is 1-2s. If we use 2s on this estimation, $5FWHM = 10s = 0.17\text{min}$. You can see the **Blank Subtracted** hits only when you click **All** in the toolbar.

Blank Hit Subtraction is performed against the “blank” sample(s). The hit(s) in any sample(s) with **Sample Type** classified as **Blank**, **DoubleBlank**, or **MatrixBlank** will automatically get subtracted from all the standard samples during the process. You can designate the “blank” sample for blank subtraction purposes by changing the **Sample Type** in the **Sample**. **No Blank Subtraction** happens if there is no “blank” sample(s). Change the sample type to turn off **Blank Subtraction**.

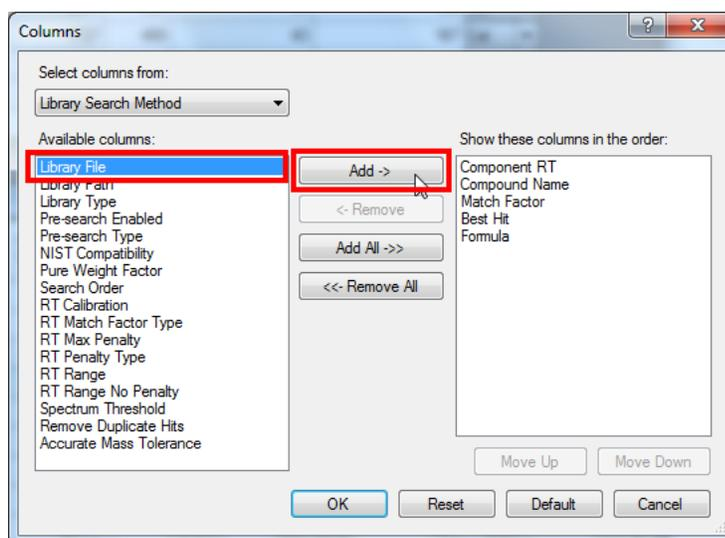
Use Show Alternate Hits to evaluate results

1. Right-click any column header in the **Components** window, and select **Add/Remove Columns**.
2. Click **Default**.



Task 2: Identify Compounds with Deconvolution

3. Select **Library File** from the **Available columns** list, and click **Add**.



4. Verify that the selected columns are moved to the **Show these columns in the order** list, and click **OK**.
5. Select the **PEST STD-10+MATRIX** sample in the **Samples** window and click **Hit** in the toolbar to view the changes in the **Components** window.

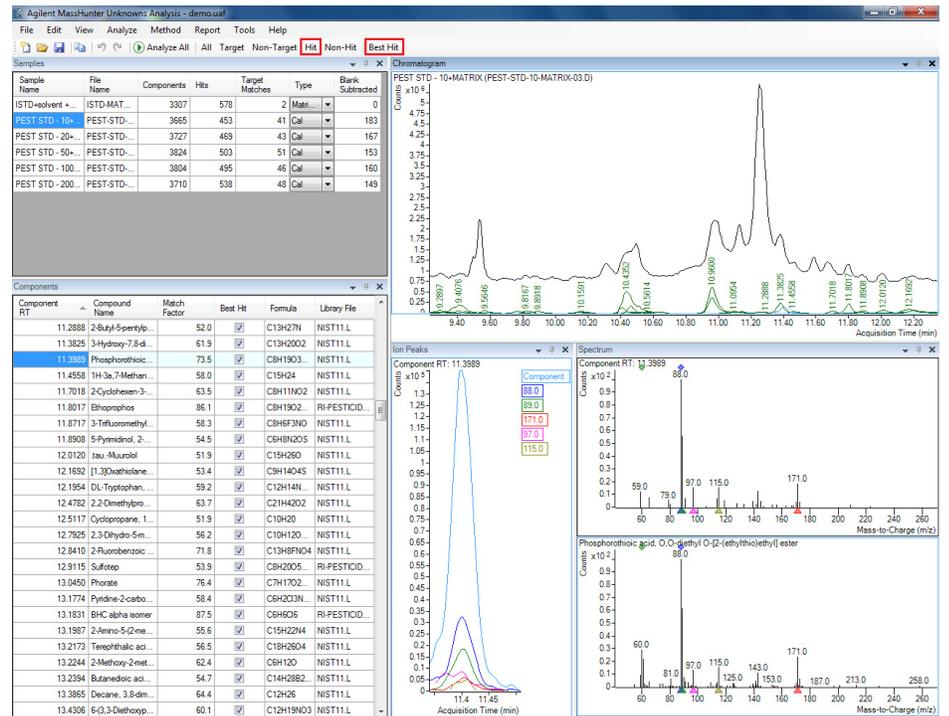
The screenshot shows the 'Agilent MassHunter Unknowns Analysis - demo.uaf' software interface. The 'Samples' table is displayed with the following data:

Sample Name	File Name	Components	Hits	Target Matches	Type
ISTD+solvent +MATRIX	ISTD-MATRI...	3307	578	2	MatrixBlank
PEST STD - 10+MATRIX	PEST-STD-1...	3665	453	41	Cal
PEST STD - 20+MATRIX	PEST-STD-2...	3727	469	43	Cal
PEST STD - 50+MATRIX	PEST-STD-5...	3824	503	51	Cal
PEST STD - 100+MATRIX	PEST-STD-1...	3804	495	46	Cal
PEST STD - 200+MATRIX	PEST-STD-2...	3710	538	48	Cal

The 'Hit' button in the toolbar is highlighted with a red box. The 'Chromatogram' window is partially visible on the right, showing a plot of 'Counts x10⁷' versus 'PEST STD - 10+MATRIX'.

Task 2: Identify Compounds with Deconvolution

Verify that the Best Hits are from different libraries.



- Right-click **Phosphorothioic acid** in the **Compounds** window and select **Show Alternate Hits**.

Component RT	Compound Name	Match
11.1417	4H-1,3-Dioxin-4-one, 2-(...	
11.2888	2-Butyl-5-pentylpyrrolidine	
11.3825	3-Hydroxy-7,8-dihydro-b...	
11.3989	Phosphorothioic acid, O...	

Add/Remove Columns...

Copy

Set Best Hits

Show Alternate Hits...

Delete Hits Del

Restore All Hits

Format Column...

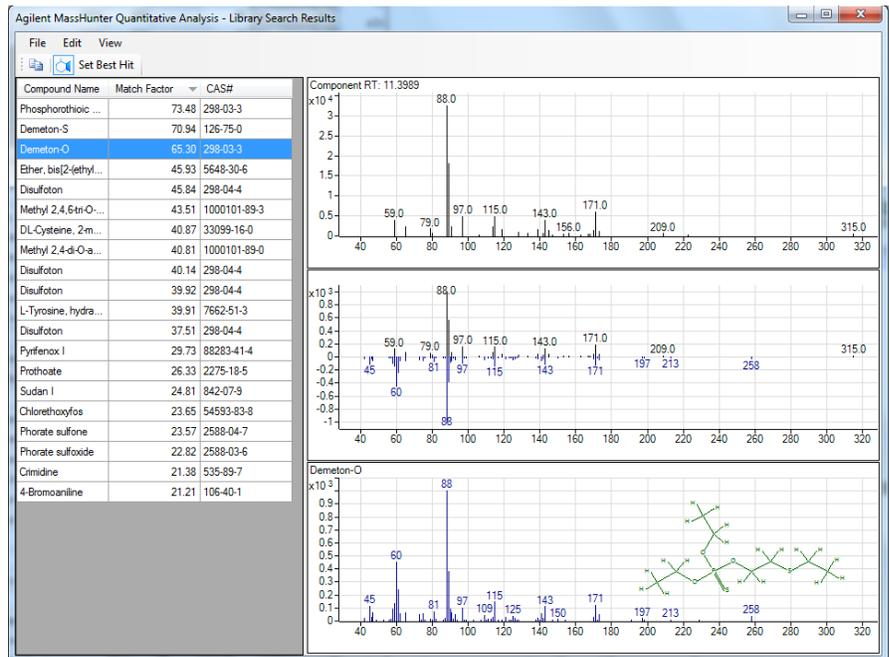
Export...

Print...

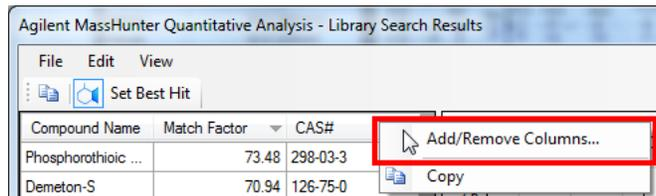
Print Preview...

Task 2: Identify Compounds with Deconvolution

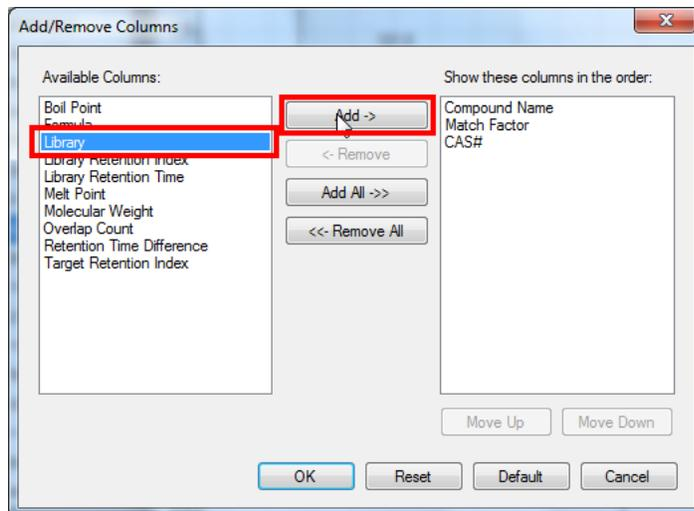
The **Library Search Results** are displayed.



- Right-click any column header in the **Library Search Results** window, and select **Add/Remove Columns**.



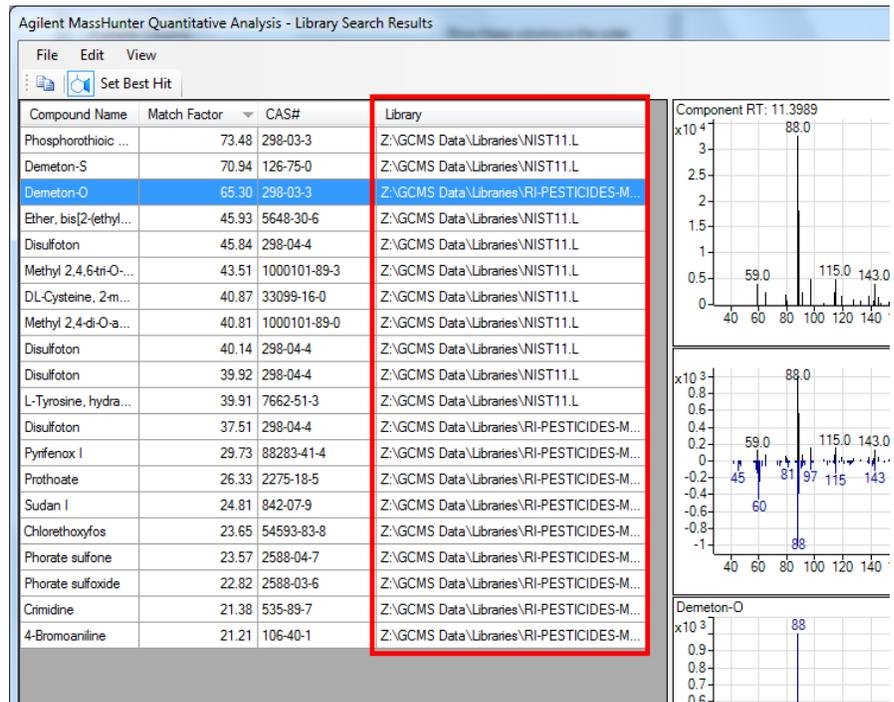
- Select **Library** from the **Available Columns** list, and click **Add**.



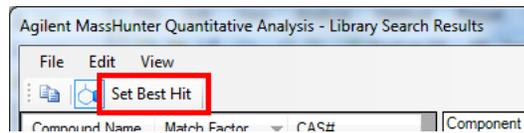
Task 2: Identify Compounds with Deconvolution

- Verify that the selected columns are moved to the **Show these columns in the order** list, and click **OK**.

The **Library** column is added to the table.



- Select **Demeton-S** and click **Set Best Hit**.



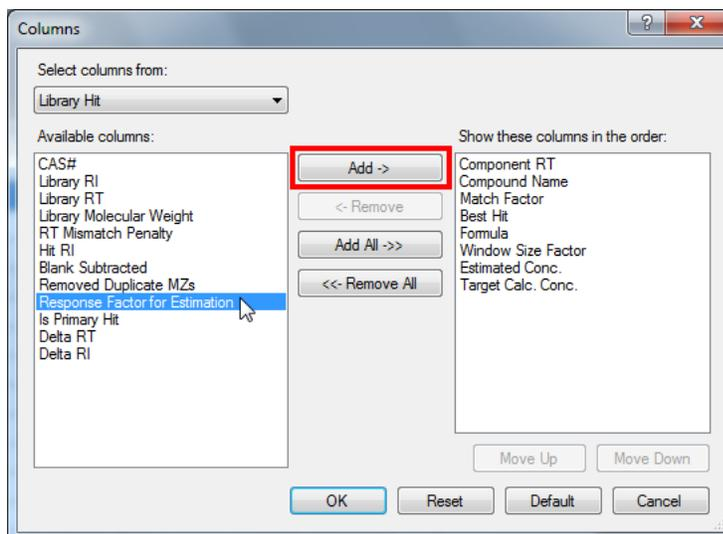
Verify that the selected compound replaced the previous compound as the current **Best Hit** in the **Component** table.

Component RT	Compound Name	Match Factor	Best Hit	Formula	Library File
11,1417	4H-1,3-Dioxin-4-one,...	50,4	<input checked="" type="checkbox"/>	C9H14O3	NIST11.L
11,2888	2-Butyl-5-pentylpyrol...	52,0	<input checked="" type="checkbox"/>	C13H27N	NIST11.L
11,3825	3-Hydroxy-7,8-dihydr...	61,9	<input checked="" type="checkbox"/>	C13H20O2	NIST11.L
11,3989	Demeton-S	70,9	<input checked="" type="checkbox"/>	C8H19O3...	NIST11.L
11,4558	1H-3a,7-Methanoaz...	58,0	<input checked="" type="checkbox"/>	C15H24	NIST11.L
11,7018	2-Cyclohexen-3-ol-1-...	63,5	<input checked="" type="checkbox"/>	C8H11NO2	NIST11.L
11,8017	Ethoprophos	86,1	<input checked="" type="checkbox"/>	C8H19O2...	RI-PESTICIDES-MOL2....
11,8717	3-Trifluoromethylben...	58,3	<input checked="" type="checkbox"/>	C8H6F3NO	NIST11.L
11,8908	5-Pyrimidinol, 2-meth...	54,5	<input checked="" type="checkbox"/>	C6H8N2OS	NIST11.L
12,0120	tau.-Muurolol	51,9	<input checked="" type="checkbox"/>	C15H26O	NIST11.L
12,1692	[1,3]Oxathiolane-4-a...	53,4	<input checked="" type="checkbox"/>	C9H14O4S	NIST11.L
12,1954	DL-Tryptophan, N-m...	59,2	<input checked="" type="checkbox"/>	C12H14N...	NIST11.L
12,4782	2,2-Dimethylpropano...	63,7	<input checked="" type="checkbox"/>	C21H42O2	NIST11.L
12,5117	Cyclopropane, 1,1,2...	51,9	<input checked="" type="checkbox"/>	C10H20	NIST11.L
12,7925	2,3-Dihydro-5-methyl...	56,2	<input checked="" type="checkbox"/>	C10H12O...	NIST11.L
12,8410	2-Fluorobenzoic aci...	71,8	<input checked="" type="checkbox"/>	C13H8FNO4	NIST11.L
12,9120	Sulfoton	71,6	<input checked="" type="checkbox"/>	C8H20O5	RI-PESTICIDES-MOL2

Task 2: Identify Compounds with Deconvolution

Review concentration estimation results

1. Right-click any column header in the **Components** window, and select **Add/Remove Columns**.
2. Select **Base Peak Deconvoluted Area, Response Factor for Estimation, Target Multiplier, Estimated Conc.,** and **Target Calc. Conc.** from the **Available columns** list, and click **Add**.
3. Verify that the selected columns are moved to the **Show these columns in the order** list, and click **OK**.
4. Select the **PEST STD-200+MATRIX** sample in the **Samples** window, and click **Target** in the toolbar to view the changes in the **Components** window.



Sample Name	File Name	Components	Hits	Target Matches	Type	Blank Subtracted
ISTD+solvent +...	ISTD-MATRI...	3307	578	2	Mat...	0
PEST STD - 10+...	PEST-STD-1...	3665	453	41	Cal	183
PEST STD - 20+...	PEST-STD-2...	3727	469	43	Cal	167
PEST STD - 50+...	PEST-STD-5...	3824	503	51	Cal	153
PEST STD - 100...	PEST-STD-1...	3804	495	46	Cal	160
PEST STD - 200...	PEST-STD-2...	3710	538	48	Cal	149

Task 2: Identify Compounds with Deconvolution

The estimated concentration results are listed in the **Estimated Conc.** column. For target compounds, you are able to compare with the Quant calculated concentrations.

Estimated Concentration is calculated using the following formula:

$$\text{Estimated Concentration} = \frac{\text{Base Peak Deconvoluted Area}}{\text{RF for Estimation}} \times \text{Multiplier}$$

Component RT	Compound Name	Match Factor	Base Peak Deconvoluted Area	Response Factor for Estimation	Target Multiplier	Estimated Conc.	Target Calc. Conc.
11.4021	Phosphorothioic aci...	94.9	2721304.4	14882.6150	1.0	182.9	144.4
11.7973	Ethoprophos	98.6	2129122.9	23030.7854	1.0	92.45	80.33
12.9031	Sulfotep	97.3	1015154.5	14085.4273	1.0	72.07	60.25
13.0452	Phorate	97.5	3420580.5	36958.8482	1.0	92.55	69.61
13.1816	BHC alpha isomer	98.8	1624103.4	30149.9823	1.0	53.87	50.11
13.7016	Pentachloroanisole	98.8	1731347.2	31680.1152	1.0	49.84	50.17
13.7651	Dimethoate	96.0	2679103.0	27597.5245	1.0	97.08	79.42
14.3218	BHC beta isomer	98.5	1014024.0	16466.2678	1.0	59.17	49.7
14.5841	Lindane	98.0	1123973.2	20441.2941	1.0	54.99	50.04
15.0133	Fonofos	97.4	2705421.8	35661.4284	1.0	75.86	60.31
15.5881	Diazinon	98.2	2028613.6	19208.4327	1.0	98.63	79.78
15.6787	Disulfoton	85.7	1203625.2	14752.9912	1.0	81.59	61.33
15.6842	BHC delta isomer	96.2	1201797.5	22136.8200	1.0	54.29	50.04
17.7267	Methyl parathion	86.9	1539303.6	7893.1866	1.0	108	71.86
17.7336	Chloropyrifos-methyl	91.9	2612062.4	25588.5789	1.0	102.1	81.29
17.9452	Heptachlor	95.7	598014.1	9867.5025	1.0	60.6	50.24

- Click **Non-Target** in the toolbar to view the estimated concentrations for Non-Targets.

The screenshot shows the Agilent MassHunter Unknowns Analysis interface. The 'Non-Target' button in the toolbar is highlighted with a red box. Below the toolbar, there is a 'Samples' table and a 'Components' table. The 'Chromatogram' window displays a plot of Counts (x10⁷) versus Acquisition Time (min) with several peaks labeled with their retention times: 11.4021, 11.7973, 12.9031, 13.0452, 13.1816, 13.7016, 13.7651, 14.3218, 14.5841, 15.0133, 15.5881, 15.6787, 15.6842, 17.7267, 17.7336, and 17.9452. The 'Components' table at the bottom shows the following data:

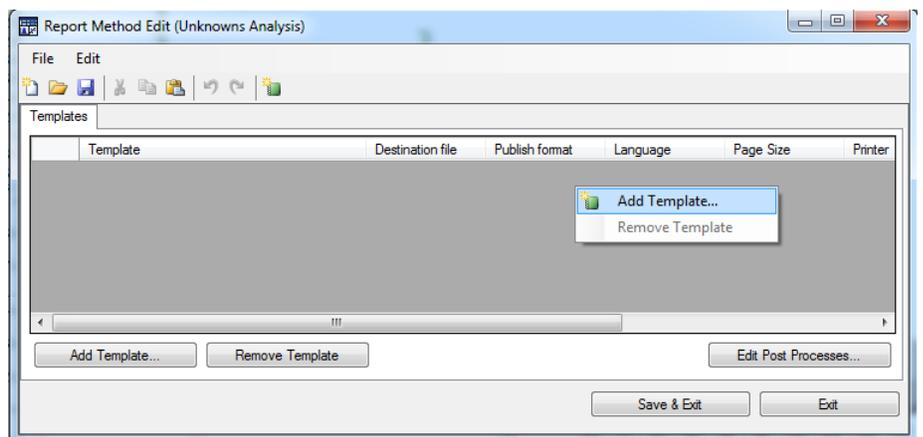
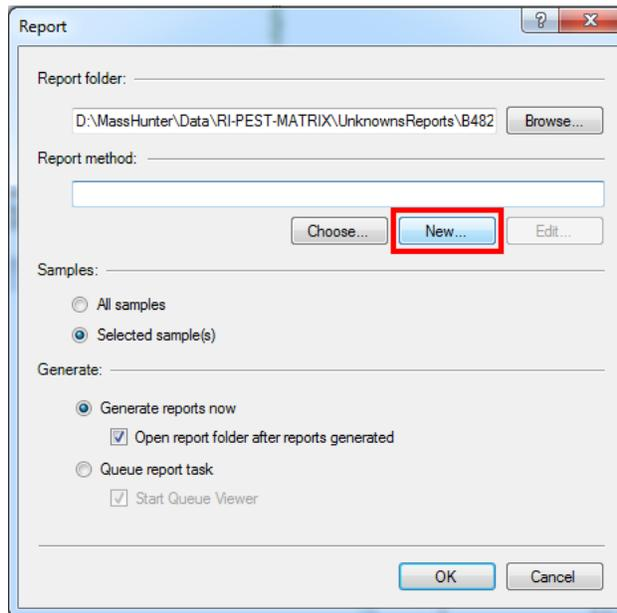
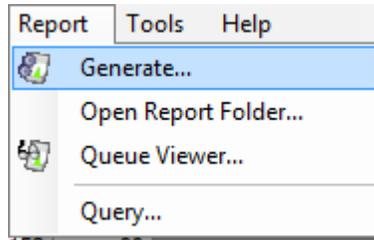
Component RT	Compound Name	Match Factor	Base Peak Deconvoluted Area	Response Factor for Estimation	Target Multiplier	Estimated Conc.	Target Calc. Conc.
37.4183	Ethanone, 1-(4-meth...	50.4	7925.2	7030154.7121		0.003202	
37.4704	Silane, diethylhepta...	77.2	73034.5	7030154.7121		0.03428	

- To save the analysis, select **File > Save Analysis**.

Task 3: Generate the Report

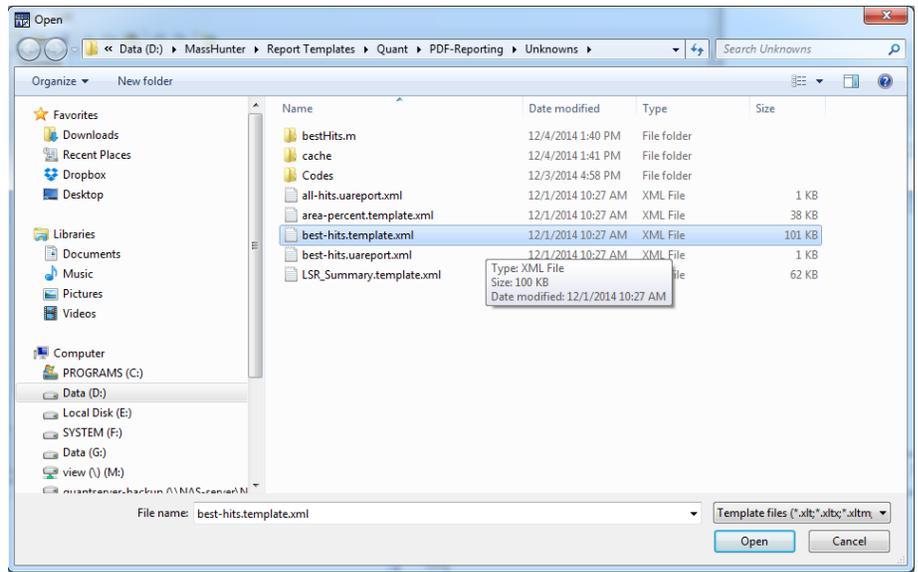
Task 3: Generate the Report

1. Select **Report > Generate**.
2. Under **Report method**, click **New**.
3. Right-click in the window and select **Add Template**.

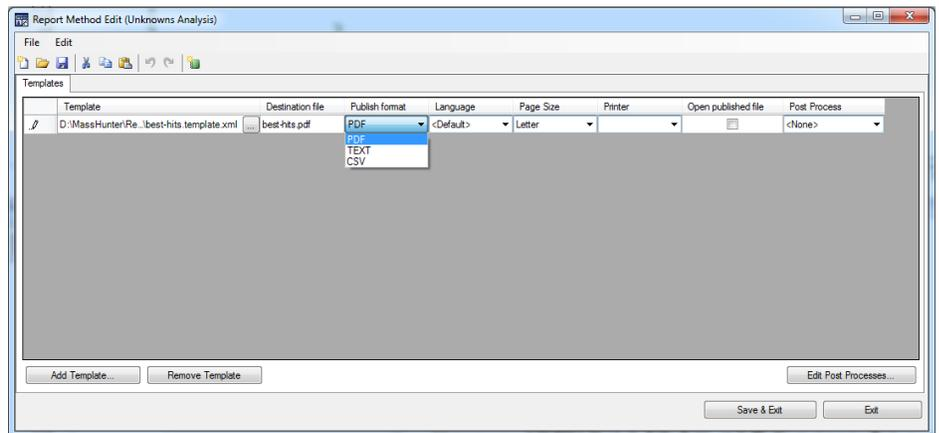


Task 3: Generate the Report

4. Navigate to **D:\MassHunter\Report Templates\Quant\PDF-Reporting\Unknowns**, select **best-hits.template.xml**, and click **Open**.

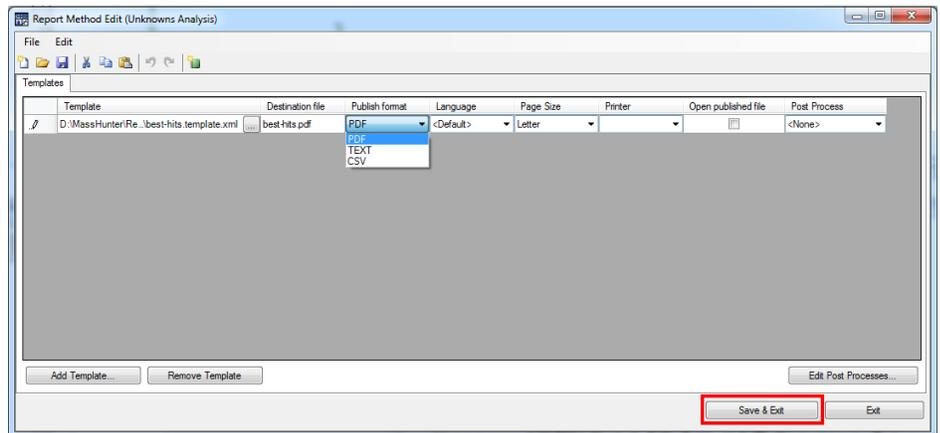


Once the template(s) is selected, you can configure the **Report Publish Format** with *PDF, TEXT, and CSV*, **Language** with *English, Chinese, Japanese, and Russian*, **Page Size**, **Printer** with *A4 and Letter*, and whether or not to **Open published file** after generating the report. The **Post Process** is also available to process the report further after finishing the report task.

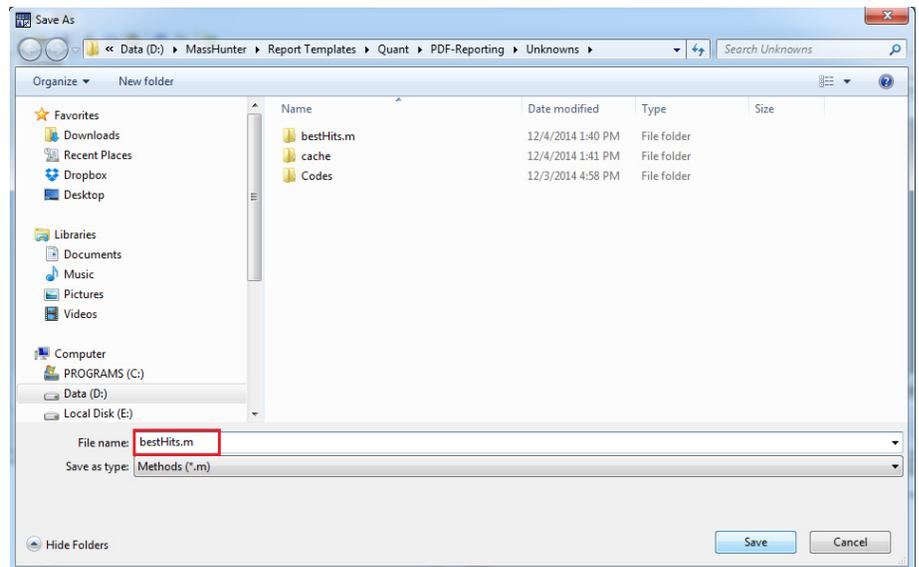


Task 3: Generate the Report

5. Click **Save & Exit** to save the Report Method in a desired location.



Report Methods have a .m extension.



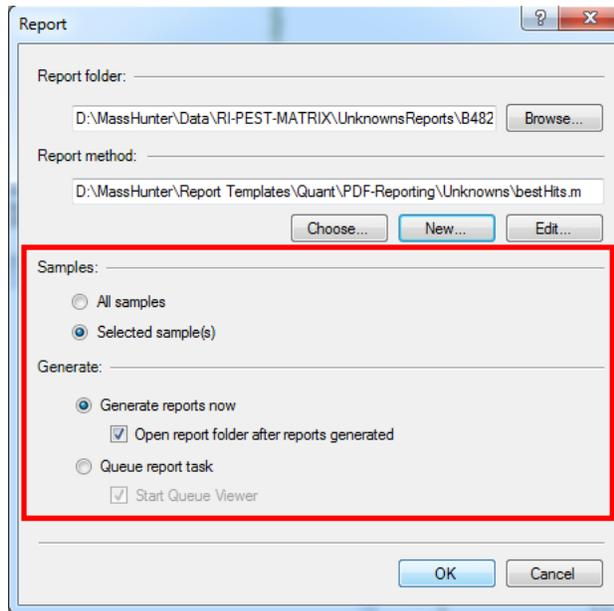
Task 3: Generate the Report

- For samples, you can generate a report for **All samples** or the **selected Sample(s)**.

For Report Generating modes, you can select **Generate reports now** or **Queue report task**.

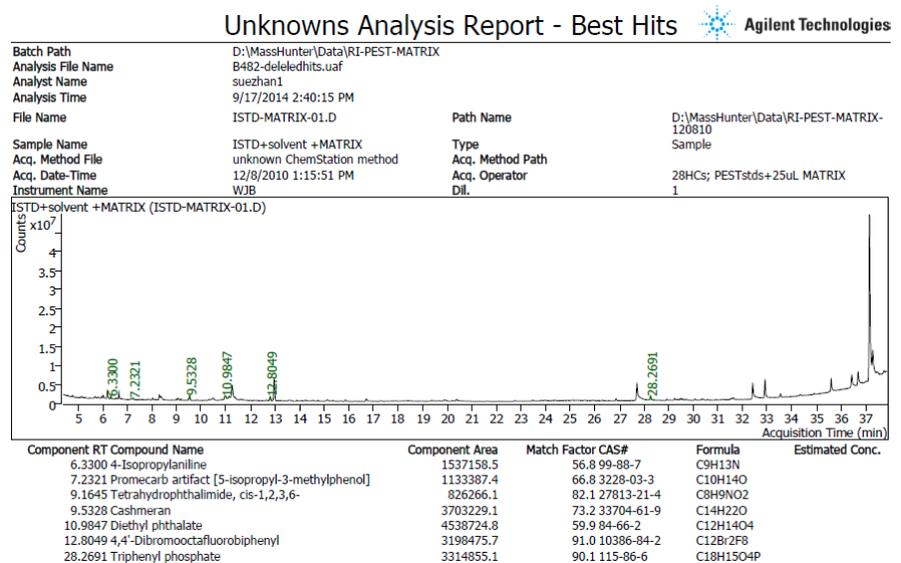
- Click **OK** to begin generating reports.

- Close the report.
- To exit the program, select **File > Exit**.



The report folder opens automatically when the report generation is complete.

Alternatively, you can select **Menu > Open Report Folder** to view the newly generated report **best-hits.pdf**. The report opens in Adobe Reader.



Task 3: Generate the Report





© Agilent Technologies, Inc.

Printed in USA, January 2015

G3335-90187