

Compound Discoverer 3.2 Metabolomics Tutorial

To familiarize yourself with the Thermo Compound Discoverer™ 3.2 application, follow the topics in this tutorial to set up a study and an analysis, process a set of example Xcalibur™ RAW files, review the result file produced by the analysis, and export the results to a Microsoft™ Excel™ spreadsheet.

Contents

- [Overview](#)
- [Start the application](#)
- [Check the computer's access to external databases](#)
- [Set up a new study and a new analysis](#)
- [Submit the analysis to the job queue](#)
- [Review the analysis results](#)
- [Modify the processing workflow and partially reprocess the analysis](#)
- [Filter the data and export the analysis results](#)

Overview

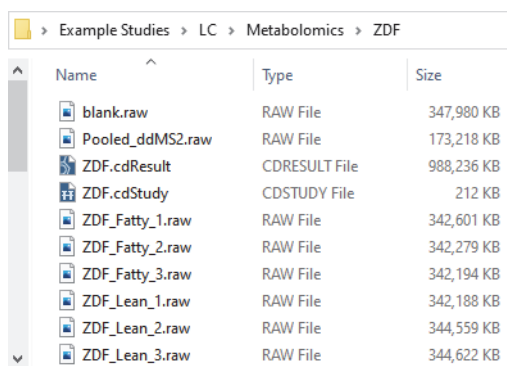
Before you open the application and begin this tutorial, review the following topics:

- [Location of the example files](#)
- [The Help system](#)
- [Tutorial summary](#)

Location of the example files

The files for this tutorial are provided on the key-shaped USB key in the software media kit. You can find these files in the following folder: Example Studies\LC\Metabolomics\ZDF.

Copy the ZDF folder to your data processing computer.



Name	Type	Size
blank.raw	RAW File	347,980 KB
Pooled_ddMS2.raw	RAW File	173,218 KB
ZDF.cdResult	CDRESULT File	988,236 KB
ZDF.cdStudy	CDSTUDY File	212 KB
ZDF_Fatty_1.raw	RAW File	342,601 KB
ZDF_Fatty_2.raw	RAW File	342,279 KB
ZDF_Fatty_3.raw	RAW File	342,194 KB
ZDF_Lean_1.raw	RAW File	342,188 KB
ZDF_Lean_2.raw	RAW File	344,559 KB
ZDF_Lean_3.raw	RAW File	344,622 KB

The Help system

The application provides Help for the application views, tabbed pages, and dialog boxes.

Do either of the following:

- From the application menu bar, choose **Help > Compound Discoverer Help**.

The Help system opens to the Welcome page.

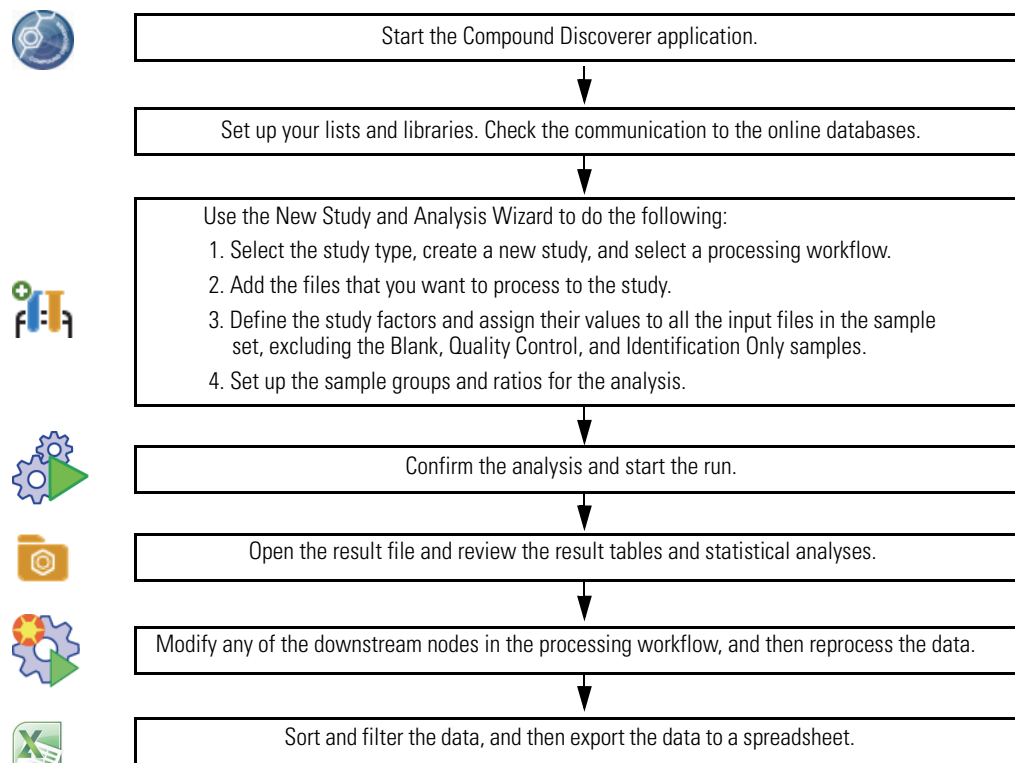
- Open a specific view, tabbed page, or dialog box. Then, on the computer keyboard, press the **F1** key.

The Help system opens to the topic for the view, page, or dialog box.



Tutorial summary

This flowchart summarizes the overall workflow for the tutorial.




Start the application

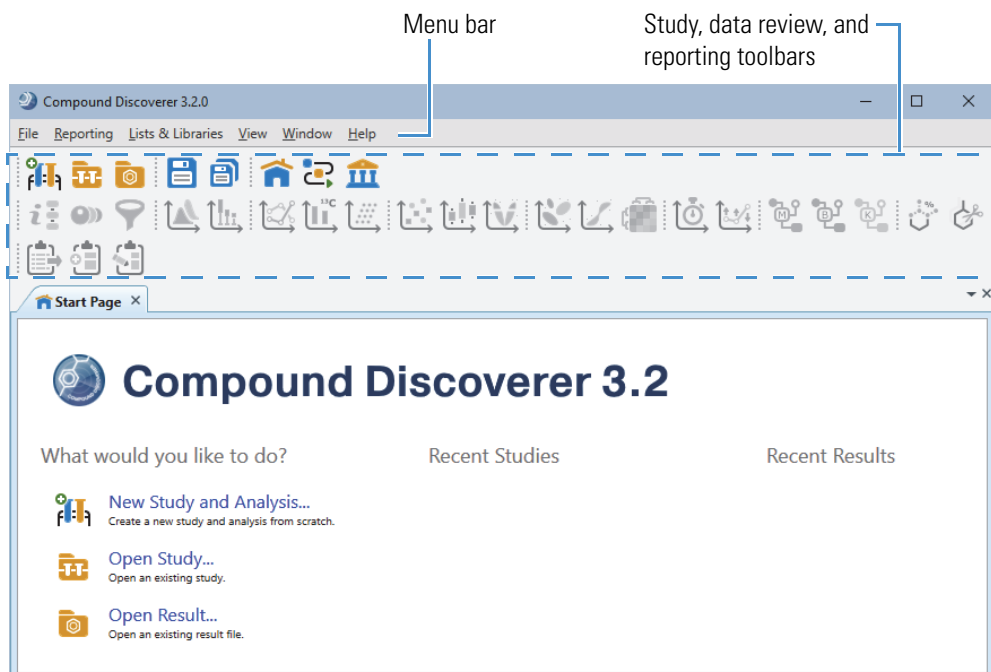
Do one of the following:

- From the taskbar, choose **Start > All Programs** (or **Programs**) > **Thermo Compound Discoverer 3.2**.

—or—

- From the computer desktop, double-click the **Compound Discoverer** icon, .

The application opens to the Start Page.



Check the computer's access to external databases

The processing workflow that you select for this tutorial uses online mass spectral databases to identify unknown compounds. To use any of the processing workflows that use online databases, such as mzCloud and ChemSpider, your processing computer must have unblocked access to these databases on the Internet.

To test the communication to the online databases that you use in this tutorial, do the following:

1. From the menu bar, choose **Help > Communication Tests**.
2. Click the **mzCloud** tab and click **Run Tests**. When the tests are complete, go to the next step.
3. Click the **ChemSpider** tab and click **Run Tests**.

If your computer has an Internet connection, but these tests fail, leave the Communication Test dialog box open and press the **F1** key to open the Help. Then, follow the instructions to troubleshoot the communication failure.

Go to the next topic to “[Set up a new study and a new analysis](#).”

Set up a new study and a new analysis

To create a new study and a new analysis, do the following:


1. [Open the New Study and Analysis Wizard](#)
2. [Select the study type, specify the directory folder, and name the new study](#)

Note The first time you set up a new study, you must specify the directory for your studies. As you create more studies, you can further organize them by creating more top-level directory folders.

3. [Select the processing workflow](#)
4. [Add the input files to the study](#)
5. [Define the experimental variables](#)
6. [Set up the sample groups and ratios](#)
7. [Customize the processing workflow](#)

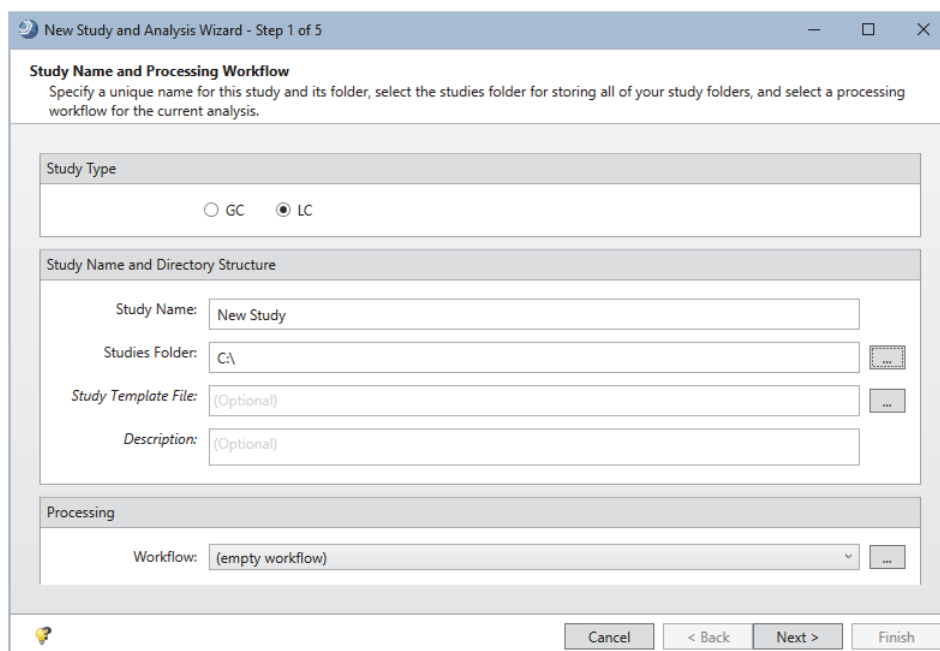
Open the New Study and Analysis Wizard

To open the New Study and Analysis Wizard, do one of the following:

- From the menu bar, choose **File > New Study and Analysis**.
- From the application toolbar, click the **Create a New Study and Analysis from Scratch** icon, .
- On the Start Page, click the **New Study and Analysis** link in the What Would You Like to Do? area.

The New Study and Analysis Wizard opens to the Study Name and Processing Workflow page.

Figure 1. Study Name and Processing Workflow page of the New Study and Analysis Wizard



Select the study type, specify the directory folder, and name the new study

Go to the next topic to “[Select the study type, specify the directory folder, and name the new study.](#)”


To select the study type, specify the directory folder, and name the new study, do the following:

1. In the Study Type area on the Study Name and Processing Workflow page of the wizard ([Figure 1](#)), select the **LC** option if it is not already selected.

The application stores this selection until you change it.

Note There are two types of studies: GC for gas chromatography-mass spectrometry data and LC for liquid chromatography-mass spectrometry data.

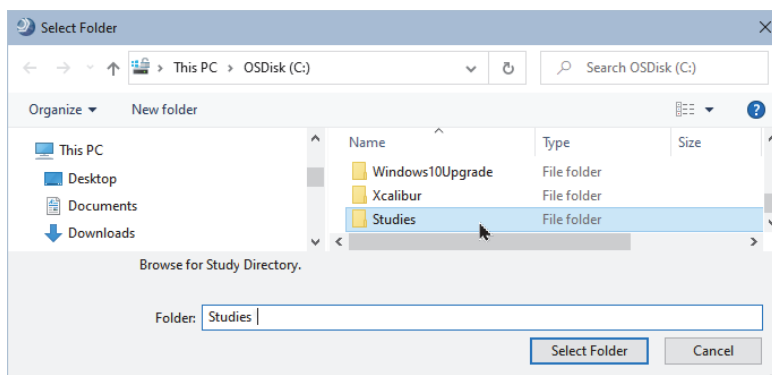
2. In the Study Name and Directory Structure area, select the studies folder as follows:

- a. Click the browse icon, , next to the Studies Folder box.

The Select Folder dialog box opens.

- b. Browse to the directory where you want to store your studies.
- c. Click **New Folder**.
- d. Name the new folder **Studies**, select it, and then click **Select Folder**.

Figure 2. Select Folder dialog box



Note The first time you create a new study, you must specify the directory (Studies Folder) where you want to store your studies. Thereafter, you can use the same studies folder or create additional studies folders.

3. In the Study Name and Directory Structure area, name the study **ZDF Example** in the Study Name box.

Note When you create a new study, the application creates a new study folder with the same name and stores the study file (.cdStudy) in the new folder and the new study folder in the specified studies folder.

Stay on this page of the wizard and go the next topic to “[Select the processing workflow.](#)”

Select the processing workflow

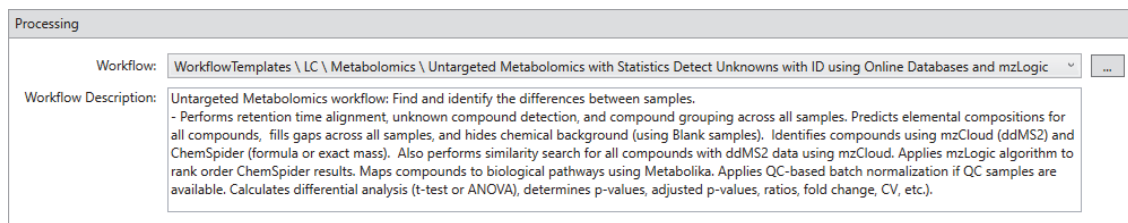
In the Compound Discoverer application, the processing method that interprets the raw data is called a processing workflow (.cdProcessingWF). The application provides defined processing workflows for several fields of study including the metabolomics field. This tutorial uses a defined processing workflow that searches the mzCloud and ChemSpider databases to identify the compounds detected in the input files (Xcalibur RAW files). This workflow also maps compounds to their biological pathways by using the local Metabolika pathway files.

To select the processing workflow, do the following:

1. In the Processing area on the Study Name and Processing Workflow page of the wizard ([Figure 1](#)), select the following processing workflow from the Workflow list:

Workflow Templates \LC\ Metabolomics\Untargeted Metabolomics with Statistics Detect Unknowns with ID Using Online Databases and mzLogic

A description of the processing workflow appears in the Workflow Description box.



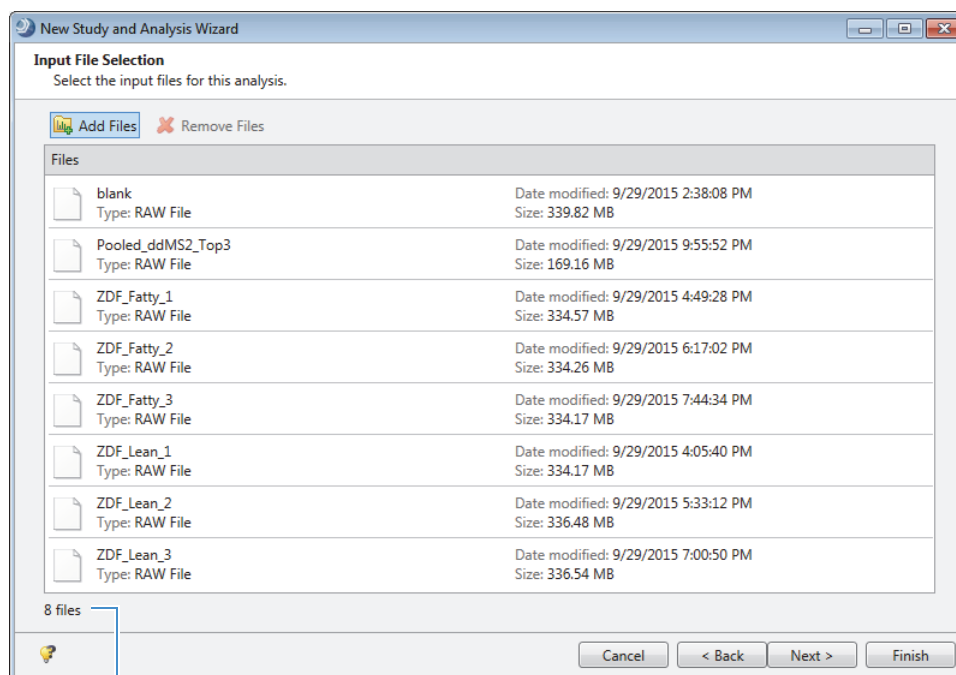
2. Read the description.
3. Click **Next** to go to the Input File Selection page of the wizard. Then, go to the next topic to [“Add the input files to the study.”](#)

Add the input files to the study

To add the input files to the study, do the following:

1. On the Input File Selection page of the wizard, click **Add Files** in the command bar. The Add Files dialog box opens.
2. Browse to the folder where you copied the Xcalibur RAW files.
3. Select all the Xcalibur RAW files in this folder and click **Open**.

The names of the selected files appear in the Files box on the Input File Selection page.

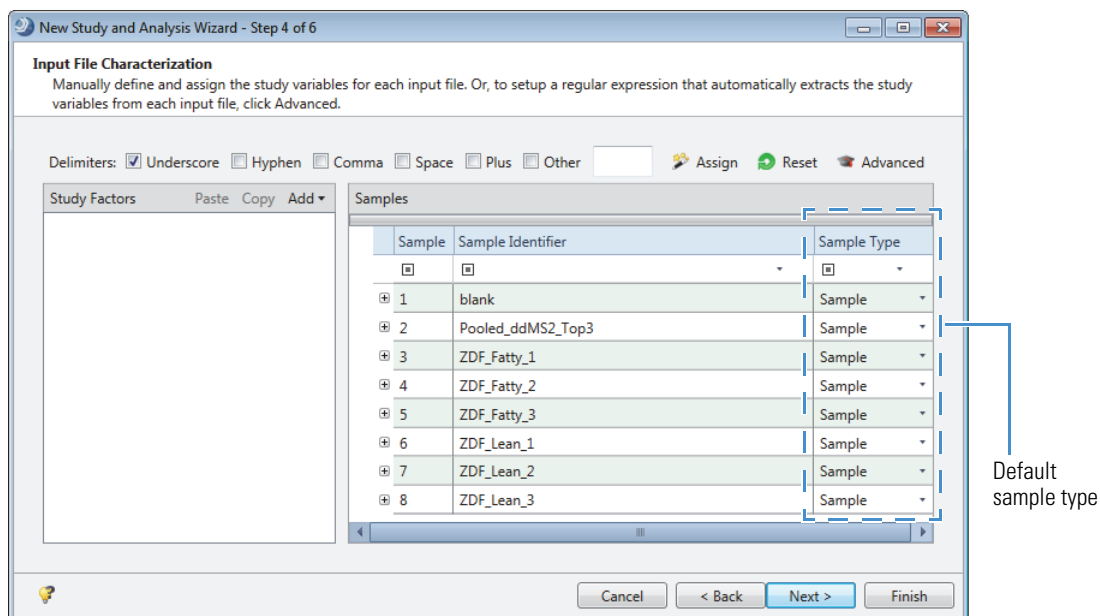


Number of files

4. Click **Next** to go to the Input File Characterization page of the wizard, and then go to the next task to [“Define the experimental variables.”](#)

Define the experimental variables

This figure shows the newly added samples in the Samples pane of the Input File Characterization page. By default, the application assigns Sample as the sample type to new samples. To group the samples according to your experimental design, you must define their study factors. In this tutorial, you are comparing the compounds in the plasma from two phenotypes of ZDF rats—Fatty and Lean.



To define and assign the study factors, do the following:

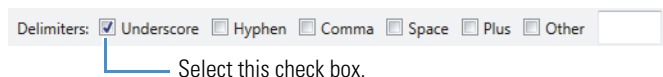
1. [Select the delimiters for parsing the file names](#)
2. [Define the phenotype as a categorical study factor](#)
3. [Assign the phenotypes and the sample types to the samples](#)

Select the delimiters for parsing the file names

In the example files, the underscore character is the delimiter that separates the study factors from the other parts of the file name. The study factor is phenotype (lean or fatty).

ZDF_Fatty_1 ZDF_Fatty_2 ZDF_Fatty_3 Blank
ZDF_Lean_1 ZDF_Lean_2 ZDF_Lean_3 Pooled_ddMS2

To select the delimiter for parsing the file names of the example files, on the Input File Characterization page, select the **Underscore** check box.



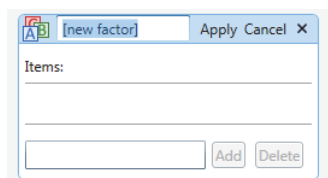
Define the phenotype as a categorical study factor

In this study, the phenotype of the Zucker diabetic fatty (ZDF) rats (fatty or lean) is a categorical study factor, as it describes an observable rather than a quantifiable feature of the group.

To define the study factor for the example study, do the following:

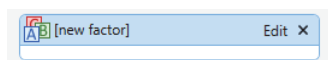
1. In the Study Factors area of the Input File Characterization page, choose **Add > Categorical Factor**.

The categorical study factor editor opens in the Study Factors area with the [new factor] box selected.

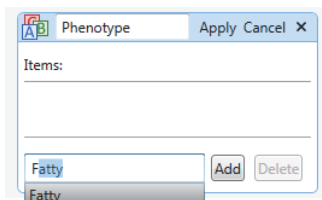


2. Type the factor name: **Phenotype**.

If the editor closes before you type Phenotype, click **Edit** to reopen it. Select [**new factor**] and type **Phenotype**.



3. To add the Fatty and Lean phenotypes to the Items list, do the following:
 - a. In the box next to the Add button, type **F**.



The application parses the file names in the Sample Identifier column and Fatty appears in the box.

- b. Click **Add**.

Fatty appears in the list.

- c. Type **L** in the box.

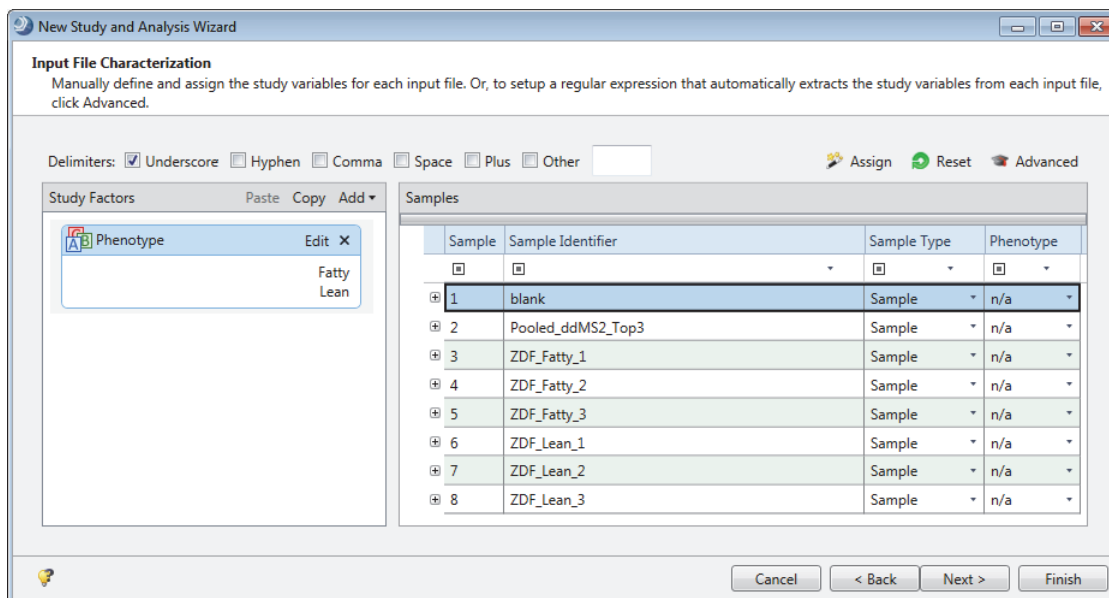
Lean appears in the box.

- d. Click **Add**.

Lean appears in the list.

4. Click **Apply**.

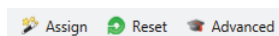
The editor closes and the Edit button replaces the Apply and Cancel buttons.



Assign the phenotypes and the sample types to the samples

To assign the study factor values and sample types to the input files, do the following:

1. In the command bar, click **Assign**.



The application does the following:

- Assigns the study factor items for the Phenotype study factor, Fatty or Lean, to the samples.
- Assigns the Blank sample type to the Blank.raw file.

This figure shows the study variable assignments. Study variables include the study factor and the sample type assignments for the samples.

Samples				
	Sample	Sample Identifier	Sample Type	Phenotype
⊞	1	blank	Blank	n/a
⊞	2	Pooled_ddMS2_Top3	Sample	n/a
⊞	3	ZDF_Fatty_1	Sample	Fatty
⊞	4	ZDF_Fatty_2	Sample	Fatty
⊞	5	ZDF_Fatty_3	Sample	Fatty
⊞	6	ZDF_Lea_n_1	Sample	Lean
⊞	7	ZDF_Lea_n_2	Sample	Lean
⊞	8	ZDF_Lea_n_3	Sample	Lean

Because the file names for the Pooled_ddMS2_Top3 sample and the Blank sample do not match the defined pattern, ZDF_categorical factor, the application assigns these samples a value of n/a for the Phenotype study factor.

IMPORTANT The assignment of n/a for a study factor is appropriate for the Blank and Identification Only sample types, as the application does not use these sample types in statistical analyses.

Note The Pooled_ddMS2_Top3 sample is a blend of the six non-blank samples. The Pooled_ddMS2_Top3.raw file acquired from this sample contains the data-dependent MS2 scans for the input file set. The full scan data for this sample was acquired at a lower resolution than that for the other samples. The Blank sample is a matrix blank and only contains the background compounds.

- For the Pooled_ddMS2_Top3 sample, select **Identification Only** in the Sample Type list.

Samples					
Error	Sampl	File	Sample Identifier	Sample Type	Phenotype
⊞	⊞	⊞	⊞	⊞	⊞
	S1	F1	blank	Blank	n/a
	S2	F2	Pooled_ddMS2	Sample	n/a
	S3	F3	ZDF_Fatty_1	Sample	Fatty
	S4	F4	ZDF_Fatty_2	Control	Fatty
	S5	F5	ZDF_Fatty_3	Blank	Fatty
	S6	F6	ZDF_Lea_n_1	Quality Control	Lean
	S7	F7	ZDF_Lea_n_2	Identification Only	Lean
	S8	F8	ZDF_Lea_n_3	Standard	Lean
				Labeled	

⌵ Show Associated File

Click **Next** to go to the Sample Groups and Ratios page of the wizard. Then, go the next topic to “[Set up the sample groups and ratios.](#)”

Set up the sample groups and ratios

Use the Sample Groups and Ratios page of the wizard to set up a differential analysis for the two phenotypes: Lean and Fatty.

To set up the sample groups and ratios, do the following:

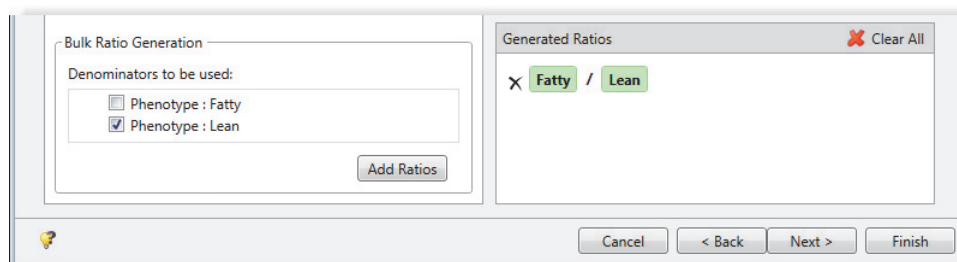
1. In the Study Variables area, select the **Phenotype** check box.

The sample groups—n/a, Fatty, and Lean—appear in the Generated Sample Groups area.

The n/a group includes only these sample types:

- Blank
- Identification Only

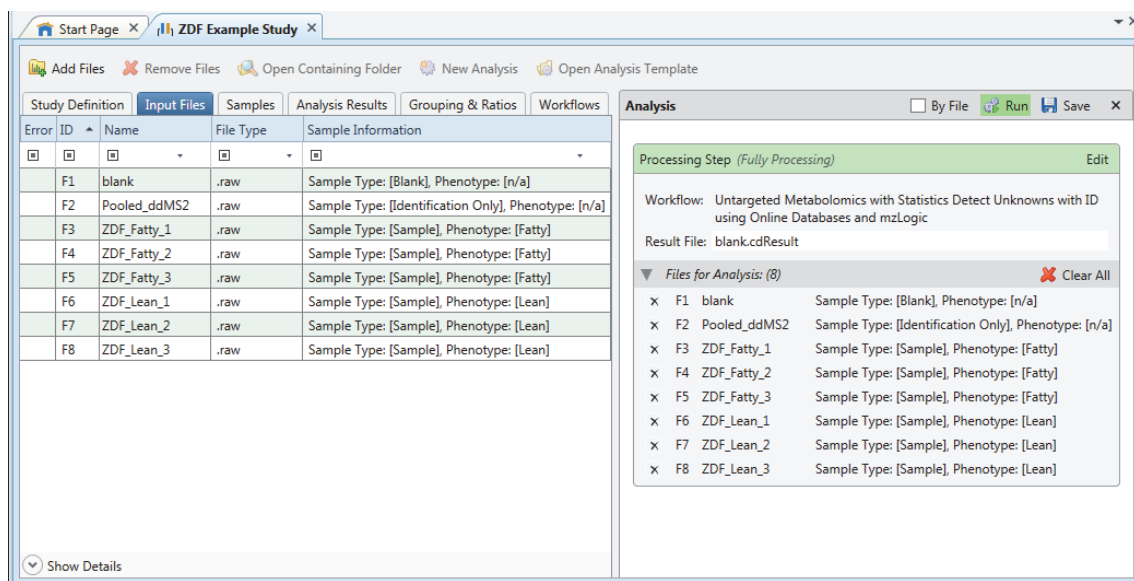
2. If an error message (❗) appears in the Generated Sample Groups area, do the following:
 - a. Return to the Input File Characterization page and verify the following sample type assignments:
 - Blank is assigned to the Blank sample.
 - Identification Only is assigned to the Pooled_ddMS2 sample.
 - b. Return to the Sample Groups and Ratios page and verify that you have corrected the sample type assignment error.
3. In the Bulk Ratio Generation area, select the **Phenotype: Lean** check box to make the Lean group the control group.
The Add Ratios button becomes available.
4. Click **Add Ratios**.
The ratio, Fatty/Lean, appears in the Generated Ratios area.
The analysis is set up to compare the two sample groups: Fatty versus Lean.



5. Click **Finish** to save the study and close the wizard.

The ZDF Example Study tab, the two analysis page tabs (Grouping & Ratios and Workflows), and the Analysis pane appear.

The Analysis pane lists the eight input files. The analysis is set up to combine the mass spectral data in these files into one result file—that is, the By File check box is clear and the file name for the result file is available for editing.



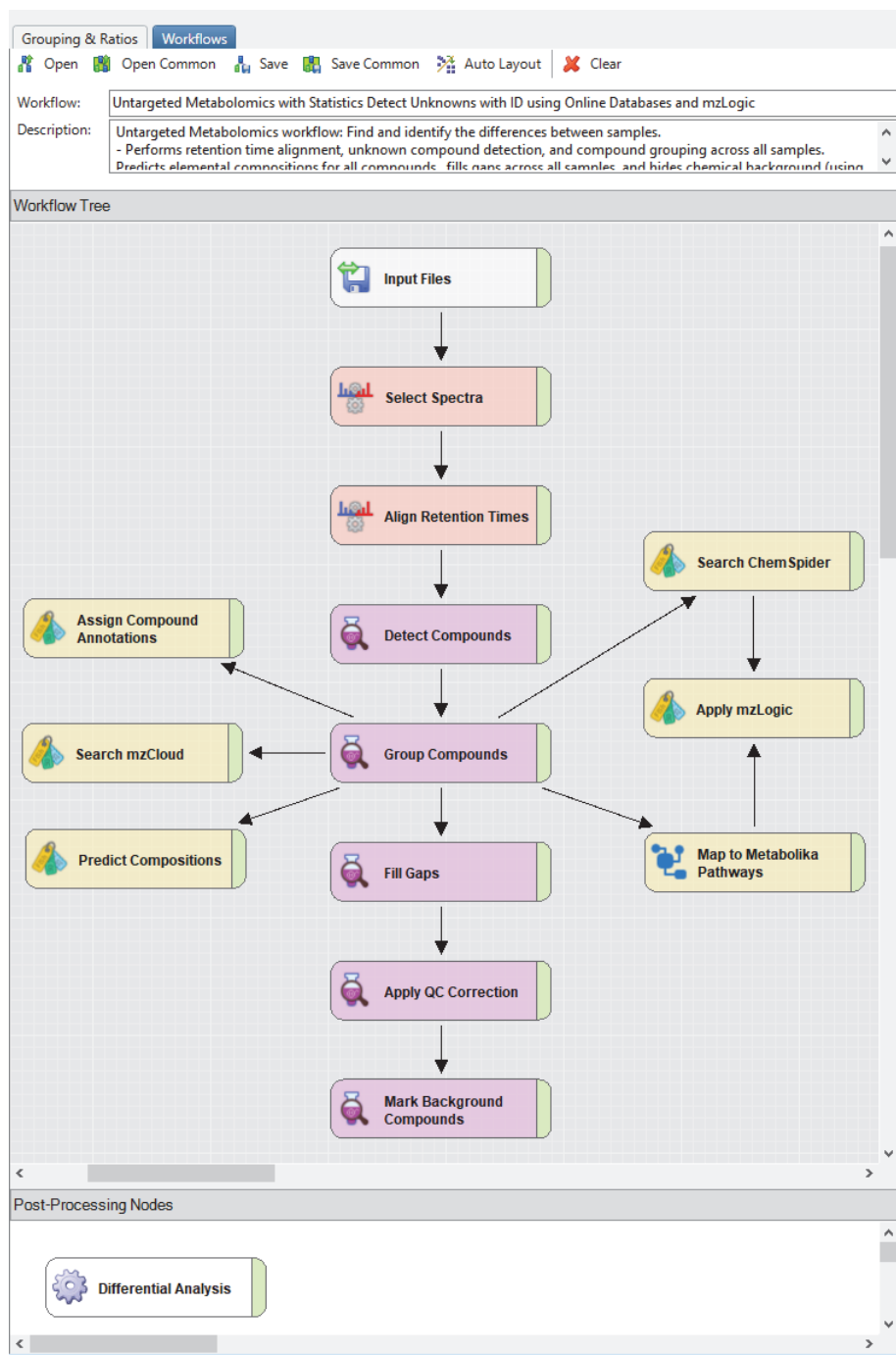
Customize the processing workflow

Before submitting the analysis to the job queue, review the processing workflow and make changes as needed.

To customize the processing workflow, do the following:

1. Click the **Workflows** tab to open the Workflows page.

Figure 3. Processing workflow for this tutorial



- Open the parameter settings for the Detect Compounds node and check the Min. Peak Intensity setting against the suggested setting for your data set (Table 1).

The minimum peak intensity setting defines the base peak intensity for the unknown compound detection. For the example ZDF data files, keep the setting of 1 000 000.

Tip To open the parameter settings for a workflow node, click the node in the Workflow Tree pane. The Parameters page for the node opens to the left.

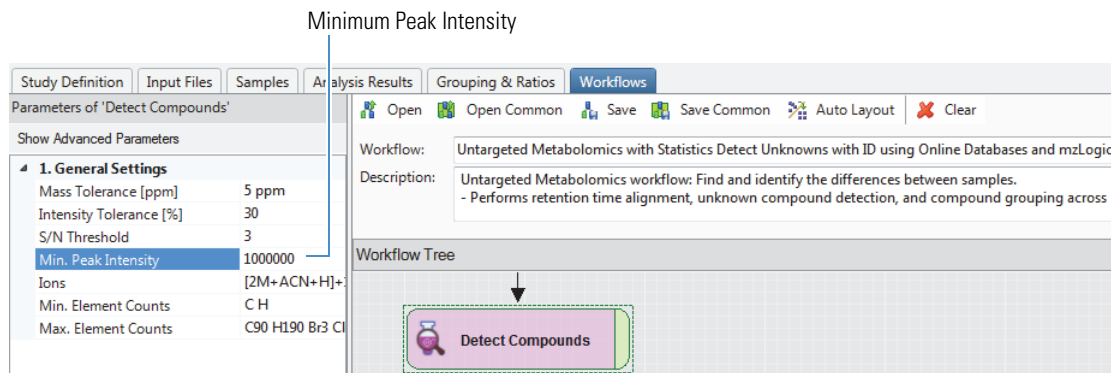


Table 1 lists the recommended range for the minimum peak intensity parameter. The optimal setting depends on the sensitivity of the mass spectrometer.

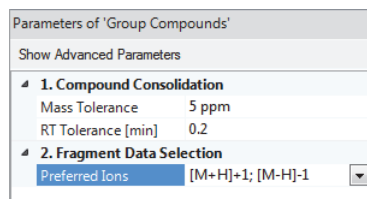
Table 1. Recommended minimum peak intensity range

Mass spectrometer	Minimum peak intensity (chromatographic peak height)
Q Exactive™, Q Exactive Plus™, Q Exactive HF	500 000 to 1 000 000
Exactive™, Exactive Plus™, Orbitrap Elite™, Orbitrap Velos Pro™	100 000 to 500 000
Orbitrap Fusion™, Orbitrap Lumos, Orbitrap ID-X	50 000 to 100 000
LTQ Orbitrap XL™, LTQ Orbitrap Velos™	25 000 to 100 000

- Open the parameter settings for the Group Compounds node.

The node is set to the default settings. For this tutorial, do not change the selection. For other analyses, select the appropriate ions from the Preferred Ions list, making sure that the selected ions include the ion selected for the Base Ions parameter in the Detect Compounds node.

For the preferred adduct ions of each detected compound, the node sends all the data-dependent MS2 scans to the search nodes and the best MS1 scan and best MS2 scan across the raw data file set to the Predict Compositions node.



IMPORTANT When setting up the node parameters, make sure that the list of preferred ions specified in the Group Compounds node includes the list of base ions specified in the Detect Compounds node.

4. Open the parameter settings for the Predict Compositions node.

The node is set to the default settings. For each detected compound, the Pattern Matching algorithm uses the best MS1 scan and the Fragments Matching algorithm uses the best MS2 scan that the analysis finds across the raw data file set.

Parameters of 'Predict Compositions'

Hide Advanced Parameters

1. Prediction Settings	
Mass Tolerance	5 ppm
Min. Element Counts	C H
Max. Element Counts	C90 H190 Br3 Cl4 N10 O18 P3 S5
Min. RDBE	0
Max. RDBE	40
Min. H/C	0.1
Max. H/C	4
Max. # Candidates	10
Max. # Internal Candidates	200
2. Pattern Matching	
Intensity Tolerance [%]	30
Intensity Threshold [%]	0.1
S/N Threshold	3
Min. Spectral Fit [%]	30
Min. Pattern Cov. [%]	90
Use Dynamic Recalibration	True
3. Fragments Matching	
Use Fragments Matching	True
Mass Tolerance	5 ppm
S/N Threshold	3

5. Open the parameter settings for the **Search mzCloud** node.

The node is set to search the entire mzCloud database, and the Identity Search is set to Cosine. For this tutorial, do not change the settings.

Parameters of 'Search mzCloud'

Hide Advanced Parameters

1. General Settings	
Compound Classes	All
Precursor Mass Tolerance	10 ppm
FT Fragment Mass Tolerance	10 ppm
IT Fragment Mass Tolerance	0.4 Da
Library	Autoprocessed; Reference
Post Processing	Recalibrated
Max. # Results	10
Annotate Matching Fragments	True
2. DDA Search	
Identity Search	Cosine
Match Activation Type	True
Match Activation Energy	Match with Tolerance
Activation Energy Tolerance	20
Apply Intensity Threshold	True
Similarity Search	None
Match Factor Threshold	60
3. DIA Search	
Use DIA Scans for Search	False
Max. Isolation Width [Da]	500
Match Activation Type	False
Match Activation Energy	Any
Activation Energy Tolerance	100
Apply Intensity Threshold	False
Match Factor Threshold	20

In the metabolomics templates, the Identity Search parameter is set to Cosine.

6. Open the parameter settings for the Search ChemSpider node.

Three out of 275 databases are selected. For this tutorial, do not change the selection.

Parameters of 'Search ChemSpider'

Hide Advanced Parameters

1. Search Settings

Database(s) BioCyc; Human Metabolome Database; KEGG

Search Mode By Formula or Mass

Mass Tolerance 5 ppm

Max. # of results per compound 100

Result Order (for Max. # of results per compound) Order By Reference Count (DESC)

Max. # of Predicted Compositions to be searched per Compound 3

2. Predicted Composition Annotation

Check All Predicted Compositions False

Show Checked Only (3/275)

Filter...

Check All Uncheck All

7. Open the parameter settings for the Map to Metabolika Pathways node.

The node is set to search all the Metabolika pathways. For this tutorial, do not change the settings.

Parameters of 'Map to Metabolika Pathways'

Show Advanced Parameters

1. Search Settings

Metabolika Pathways \\(3R)-linalool biosynthesis.metabolika

Search Mode By Formula or Mass

2. By Mass Search Settings

Mass Tolerance 5 ppm

3. By Formula Search Settings

Max. # of Predicted Compositions to be searched per Compound 3

4. Display Settings

Max. # Pathways in 'Pathways' column 20

Submit the
analysis to
the job queue

When the Run command in the Analysis pane becomes available, you can start the processing run.

To rename the file name of the result file and submit the input files for processing together into one result file, do the following:

1. To create one result file for the input file set, leave the **By File** check box clear.
By default, the application uses the name of the first input file as the result file name.
2. In the Result File box, change file name for the result file to **Fatty vs Lean**.

Analysis

By File Run Save X

Processing Step (Fully Processing) Edit

Workflow: Untargeted Metabolomics with Statistics Detect Unknowns with ID using Online Databases and mzLogic

Result File: Fatty vs Lean.cdResult

Run command


Result file name

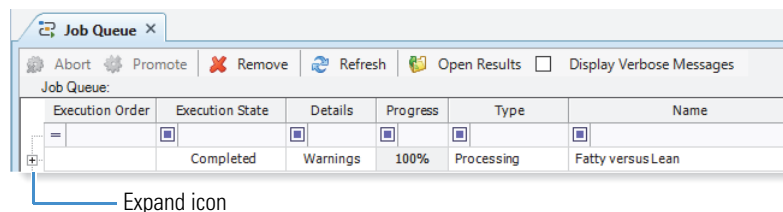
3. To submit the analysis to the job queue, click **Run**.

Tip The processing workflow for this tutorial includes the Differential Analysis node. If no groups or ratios are set up on the Grouping & Ratios page of the analysis, a warning prompt appears. If you are not interested in grouping or comparing the samples, you can ignore this warning.

4. If a warning prompt appears while you are working with this tutorial, do the following:
 - a. Click **Abort**.
 - b. Open the **Grouping & Ratios** page.
 - c. In the Study Variables area, select the **Phenotype** check box if it is not already selected.
 - d. In the Bulk Ratio Generation area, select the **Lean** check box if it is not already selected and click **Add Ratios**.
 - e. To restart the analysis, click **Run** in the Analysis pane.

The Job Queue page opens.

- To view the processing messages, click the expand icon, , to the left of the job row.



Note During the run, the Map to Metabolika node and the Apply QC Correction node generate warning messages, which you can ignore.

Warning messages have a yellow background.

(34): Map to Metabolika Pa...	Warning	8 pathway compounds with invalid elemental composition not persisted.
(39) Apply QC Correction	Warning	Cannot find any QC sample. QC correction will be skipped.

Review the analysis results

To review the analysis results, follow these topics in order or jump to the topics of interest:

- [Open the result file](#)
- [Review the default result page layout](#)
- [Modify the layout of the result tables or views](#)
- [Review the chromatograms for each compound](#)
- [Review the full scan MS1 spectra](#)
- [Review the fragmentation spectra](#)
- [Apply the Statistics layout](#)
- [View a trend chart for a single compound or a set of trend lines for multiple compounds](#)
- [Visualize the detected compounds as a volcano plot](#)
- [View the principal component analysis](#)
- [Find a set of discriminating compounds](#)
- [Run a hierarchical clustering analysis](#)
- [View the Metabolika pathways for a compound](#)
- [Review the results of the mzLogic analysis](#)

Open the result file

You can open a result file from multiple locations: the Job Queue page, the Analysis Results page of a study, the Compound Discoverer Start Page, or the menu bar.

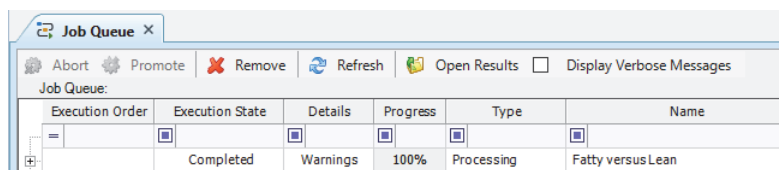
Tip If you did not reprocess the example data set, do the following:

1. Open the result file provided on the Compound Discoverer 3.2 USB key:

Example Studies\Metabolomics Study\ZDF Study\ZDF.cdResult

2. From the menu bar, choose **Window > Reset Layout**.
3. Sort the Compounds table in descending order by the Area (Max) column.

To open the result file when the run is completed, double-click the run on the Job Queue page.



Tip If the Job Queue page is closed, from the menu bar, choose **View > Job Queue** to open it.

Review the default result page layout

The factory default layout for a result file includes the following items:

- A page tab with the result file name.
- A Chromatograms view on the top left that is populated with XIC traces (across the input file set) for the compound in the first row of the active table. The view automatically zooms in to the start and end points of the chromatographic peak for the compound, and the integrated peak areas are shaded.
- A Mass Spectrum view on the top right that is populated with the MS1 scan (for a preferred ion) that is closest to the chromatographic peak apex for the selected compound across the input files. The spectrum tree to the left includes the MS1 scans and the fragmentation scans for the preferred ions that were acquired within the following time range:
 - The chromatographic peak apex for the selected compound \pm peak width at half maximum (FWHM)
- or–
- The Start and end points of the chromatographic peak, as determined by the peak detection algorithm

Note If the data set does not include data-dependent MS2 scans within the retention time window but does include DIA scans within this window, the spectrum tree includes the DIA scans. Your Thermo Scientific data-visualization application displays DIA or AIF in the scan headers of data-independent scans. The Compound Discoverer application displays DIA for both of these scan types.

- A set of tabbed main tables below the two graphical views. For a metabolomics analysis, the Compounds table is the active table and is sorted by the Area (Max.) column—that is, the first row displays the compound with the largest chromatographic peak area (found in one of the input files).

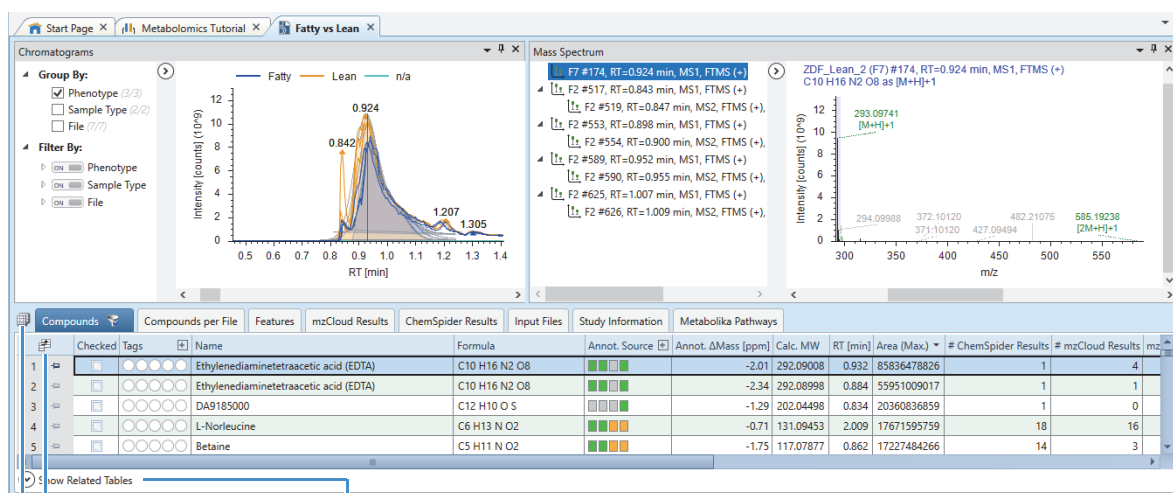
Note Because the selected processing workflow includes the Mark Background Compounds node, the Compounds tab has a filter icon with a check mark (☑). The compounds that the analysis identified as background compounds are marked as background compounds in both the blank and non-blank samples and are hidden from the table.

- A collapsed area for the related tables below the main tables.

Figure 4 shows the factory default layout for the Fatty vs Lean.cdResult file (or the ZDF.cdResult file).

Note In Compound Discoverer 3.1 or later, the chromatographic peak area for compounds and features is reported in counts \times seconds. The chromatographic peak area for a compound in each input file is the combined peak area for all its detected preferred ions (features).

Figure 4. Default result file layout



Opens the pane of related tables.

Field Chooser icon

Select Table Visibility icon

Table 2 describes the main result tables that the selected processing workflow produces.

Table 2. Main tables and some of the related tables for the selected processing workflow

Result table	Description
Visible main (top-level) tables	
Compounds	Lists all the compounds that the analysis detected, grouped by their molecular weight and retention time (MW×RT) dimensions across all the input files. The RT of a compound is the weighted average (by area) of the compound across all the input files (see the related Compounds per File table)
Compounds per File	Lists all the compounds that the analysis detected across all the input files on a per file basis. Does not list compounds that the Fill Gaps node detected by filling a full gap.
Features	Lists all the features (ions with the same mass-to-charge values and retention time) that the analysis detected across all the input files on a per file basis.
mzCloud Results	Lists the mzCloud search results across all the input files.
ChemSpider Results	Lists the ChemSpider search results across all the input files.
Input Files	Describes the input files that the application processed to create the result file.
Study Information	Displays the file name, sample type, and study factor values for each input file. Also displays the ratios that include the input file.
Metabolika Pathways	Lists the Metabolika Pathways that contain at least one of the detected compounds across the input files.
Visible tables related to the selected compound in the Compounds table	
Structure Proposals	Displays your structure proposals for the selected compound. Initially, this table is empty.
Compound per File	Displays information about the selected compound on a per file basis.
Predicted Compositions	Displays the predicted compositions for the selected compound. The predicted compositions are based on the molecular weight of the neutral compound.
Metabolika Results	Displays information about the matched and partially matched structures for the selected compound.
mzCloud Results	Displays the mzCloud results for the selected compound.
ChemSpider Results	Displays the ChemSpider results for the selected compound.
Metabolika Pathways	Lists the pathways that include the selected compound.
Visible table related to the selected feature in the Features table	
Chromatogram Peaks	Describes the chromatographic peak for the selected feature.
Visible table related to the selected file in the Input Files table	
File Alignments	Describes the alignment for the input file selected in the Input File table.
Hidden tables	
Adducts (main table)	Lists the adducts in the Adducts library.
Filled Gaps (related table)	Provides information about the imputed chromatographic peak areas.


Modify the layout of the result tables or views

Select the Visibility of the Result Tables

To learn how to modify the layout of the result tables, do the following:

- [Select the Visibility of the Result Tables](#)
- [Show the related tables](#)
- [Show or hide table columns](#)
- [Display the input file names for the subcolumns](#)
- [Sort the rows in a result table](#)
- [Pin a row to the top of the result table](#)
- [Freeze a table column](#)
- [Drag a view to another location or monitor](#)

The default layout does not include all the results tables that the selected processing workflow produces. By default, the Filled Gaps related table produced by the Fill Gaps node is a hidden table in the first set of related tables.

1. Click the **Select Table Visibility** icon, , to the left of the result table tabs.

The Select Visible Tables dialog box opens.



2. To display a table, select its check box. To hide a table, clear its check box.
For this tutorial, do not change the selections.
3. Click **OK**.

Show the related tables

The main tables provide information about each compound across the input file set. The tables related to a main table provide details about the item selected in the main table. For example, the Compounds per File table for a compound in the main Compounds table provides information about the compound in each input file where it is detected.


To view the tables related to the Compounds table, click **Show Related Tables** at the bottom of the page.

Tip To view the information for a specific compound in the main Compounds table, do the following:

1. Select the compound in the main table.
2. Click **Show Related Tables**.
3. In the related tables pane, click the tab of interest.

Show or hide table columns

To show or hide columns in the Compounds table, do the following:

1. Click the **Field Chooser** icon, , to open the Field Chooser box.

In the Compounds table, the following columns are hidden:

- #Adducts
- #Metabolika Pathways
- #Similarity Results
- Annot. Δ Mass [ppm]
- Annotation MW
- Area (per input file)
- Background
- FISH Coverage
- Gap Fill Status
- Gap Status
- Metabolika Pathways
- MS Depth
- RT Tolerance [min]
- Structure

2. Do one of the following:


- To display a column, select its check box.


For example, do display the Area column in the Compounds table, select the Area check box.

The Area column appears in the Compounds table to the left of the Group Areas column.

Area							Group Areas	
5.99e10	4.78e10	5.39e10	7.55e10	8.59e10	7.01e10	1.40e5	5.39e10	7.55e10
1.23e10	2.04e10	1.86e10	4.37e9	2.54e9	5.40e9	6.83e4	1.86e10	4.37e9
1.25e10	1.77e10	1.67e10	9.63e9	8.73e9	1.08e10	7.31e5	1.67e10	9.63e9

- To hide a column, clear its check box.

3. To close the Field Chooser box, click its close icon, .


Field Chooser 

<input type="checkbox"/>	# Adducts
<input checked="" type="checkbox"/>	# ChemSpider Results
<input type="checkbox"/>	# Metabolika Pathways
<input checked="" type="checkbox"/>	# mzCloud Results
<input type="checkbox"/>	# Similarity Results
<input checked="" type="checkbox"/>	Adj. P-value
<input checked="" type="checkbox"/>	Annot. Source
<input type="checkbox"/>	Annot. Δ Mass [Da]
<input checked="" type="checkbox"/>	Annot. Δ Mass [ppm]
<input type="checkbox"/>	Annotation MW
<input type="checkbox"/>	Area
<input checked="" type="checkbox"/>	Area (Max.)
<input type="checkbox"/>	Background
<input checked="" type="checkbox"/>	Calc. MW
<input checked="" type="checkbox"/>	Checked
<input type="checkbox"/>	FISH Coverage
<input checked="" type="checkbox"/>	Formula
<input type="checkbox"/>	Gap Fill Status
<input type="checkbox"/>	Gap Status
<input checked="" type="checkbox"/>	Group Areas
<input checked="" type="checkbox"/>	Group CV [%]
<input checked="" type="checkbox"/>	Log2 Fold Change
<input type="checkbox"/>	Metabolika Pathways
<input type="checkbox"/>	MS Depth
<input checked="" type="checkbox"/>	MS2
<input checked="" type="checkbox"/>	mzCloud Best Match
<input checked="" type="checkbox"/>	mzCloud Best Match Confidence
<input type="checkbox"/>	mzCloud Library Matches
<input checked="" type="checkbox"/>	Name
<input checked="" type="checkbox"/>	P-value
<input checked="" type="checkbox"/>	Ratio
<input checked="" type="checkbox"/>	RT [min]
<input type="checkbox"/>	RT Tolerance [min]
<input type="checkbox"/>	Structure
<input checked="" type="checkbox"/>	Tags

Display the input file names for the subcolumns

In the Compounds table, the following columns contain a subcolumn for each input file in the processed data set: Metabolika Pathways, Areas (hidden by default), Group Area, Group CV [%], Ratio, Log2 Fold Change, P-Value, and Adj. P-Value.

To display the headings for the subcolumns in a table column, do the following:

1. Click the expand icon, , to the right of the column name.

The names of the input files appear vertically above the individual area columns. The name of the first input file, blank.raw (F1), appears in bold font and has an asterisk indicating that it is the selected area column.

Area							Group Areas	
<div> <div>Selected column</div> <div> <div>ZDF_Fatty_1.raw (F3)*</div> <div>ZDF_Fatty_2.raw (F4)</div> <div>ZDF_Fatty_3.raw (F5)</div> <div>ZDF_Lean_1.raw (F6)</div> <div>ZDF_Lean_2.raw (F7)</div> <div>ZDF_Lean_3.raw (F8)</div> <div>blank.raw (F1)</div> </div> <div>Collapse icon</div> </div>								
5.99e10	4.78e10	5.39e10	7.55e10	8.59e10	7.01e10	1.40e5	5.39e10	7.55e10
1.23e10	2.04e10	1.86e10	4.37e9	2.54e9	5.40e9	6.83e4	1.86e10	4.37e9
1.25e10	1.77e10	1.67e10	9.63e9	8.73e9	1.08e10	7.31e5	1.67e10	9.63e9

Sort the rows in a result table


2. Click the collapse icon, , to hide the sample names.
3. To hide the Area column for the remaining portion of this tutorial, open the Field Chooser box and clear the **Area** check box.

Table 3 describes how to sort the rows in a result table.

Note The application treats formulas the same as text strings and sorts them by the order of the characters in the formula string, not by the actual number of elements in the formula.

Table 3. Sorting the rows in a result table

To do this	Do the following
Sort a result table by a column with numeric or text information.	Click a column header once or twice to sort the rows in ascending order (▲) or descending order (▼), based on the contents of the column.
Sort a result table by a column that contains a distribution map.	<ol style="list-style-type: none">1. Click the expand icon to display the vertical headings of the subordinate columns.2. Select the heading of the subordinate column that you want to sort by. The selected subordinate column heading appears in bold text.3. Click the column heading to sort the table rows.
Sort a result table by multiple columns.	<ol style="list-style-type: none">1. Click the column header of the primary sort column once or twice to sort the rows in ascending order (▲) or descending order (▼), based on the contents of the column.2. Hold down the CTRL key and click the column header of the secondary sort column once or twice to set the sort order.

Pin a row to the top of the result table



To freeze a table row to the top of the result table, click the pin icon to the right of the row number (unpinned, , or pinned, )

Figure 6 on page 21 shows row 8 (creatine) pinned to the top of the Compounds result table.

Freeze a table column

To freeze a table column to the leftmost position, do the following:

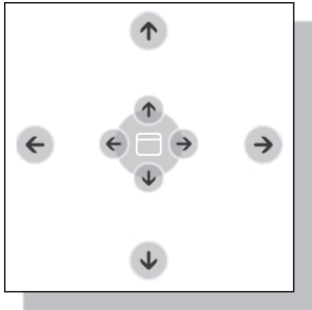
1. Right-click the result table and choose **Enable Column Fixing**.
2. Click the pin icon to the right of the column heading.

Use the mouse to move a view or tabbed document to another position as follows:

1. Drag the view by its title bar or the document by its tab.

As you drag a view or a tabbed page by its title bar, a guide tool appears. The guide tool consists of four directional arrows (inner arrows) that are arranged in a diamond pattern around a central circle. In addition to the guide tool, a directional arrow (outer arrows) appears in the middle of each of the four window edges.

Figure 5. Guide tool



2. After the guide tool appears, align the pointer with the appropriate directional arrow, and then release the mouse button.

You can also drag views to another monitor.

Open or close a view

Review the chromatograms for each compound

Modify the display options

To open or close a view, do the following:

- To open a view, choose **View > Specific view** from the menu bar or click the toolbar icon for the view.
- To close a view, click the close icon in the upper right corner of the view.

With the factory default layout, the Chromatograms view displays the overlaid traces for the compound that you select in the Compounds table, and the integrated area of each chromatographic peak is shaded.

Follow these topics to modify the display options and view specific traces in the Chromatograms view:

- [Modify the display options](#)
- [Filter the chromatographic traces](#)

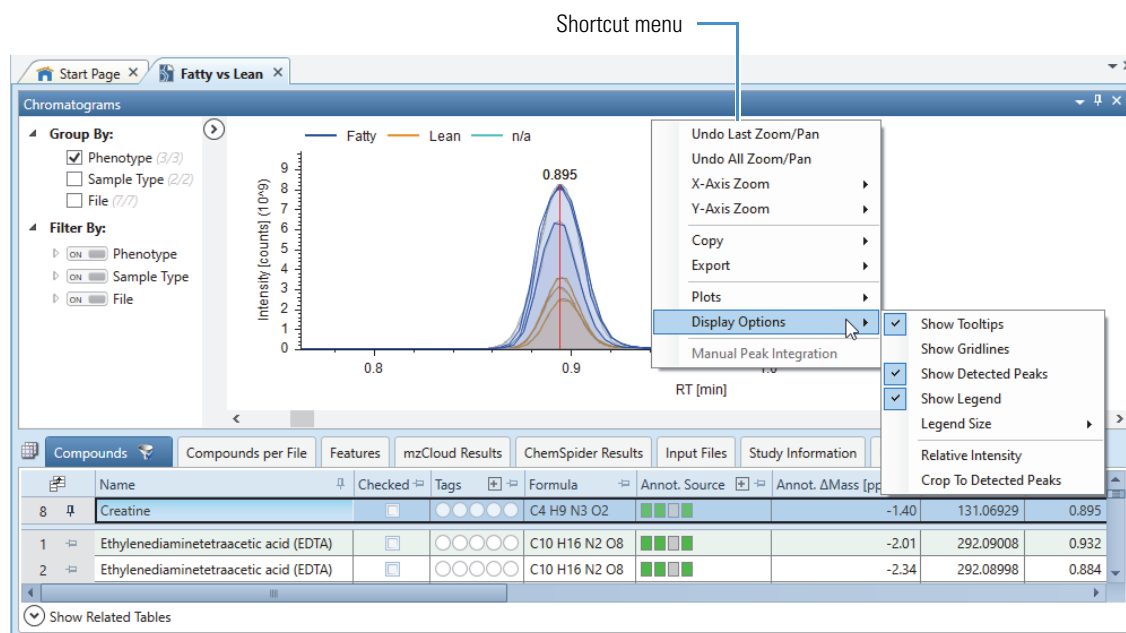
To view the chromatographic traces without the shading that can mask underlying peaks, do the following:

1. Sort the Compounds table by Area (Max.) in descending order.
2. Select **row 8** to display the overlaid XIC traces for creatine.

[Figure 6](#) displays the overlaid traces for the input files that contain the selected compound as well as the shortcut menu for the Chromatograms view. By default, the display options for the Chromatograms view are set to Show Tooltips, Show Detected Peaks, and Show Legend.

In [Figure 6](#), row 8 (creatine) is pinned the top of the Compounds result table, and the Mass Spectrum view is closed.

Figure 6. Shaded and overlaid chromatographic peaks for creatine—the compound selected in the Compounds table

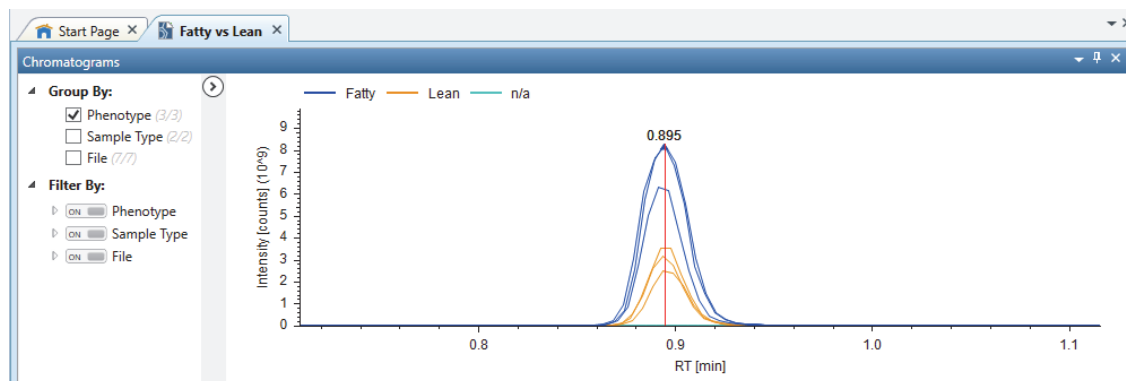


3. Right-click the Chromatograms view and choose **Display Options > Show Detected Peaks**.

Figure 7 shows the overlaid traces for all the samples.

The application displays the three sample groups in different colors. The legend shows the display colors for the groups—blue for Fatty, and orange for Lean, and turquoise for N/A (the blank sample).

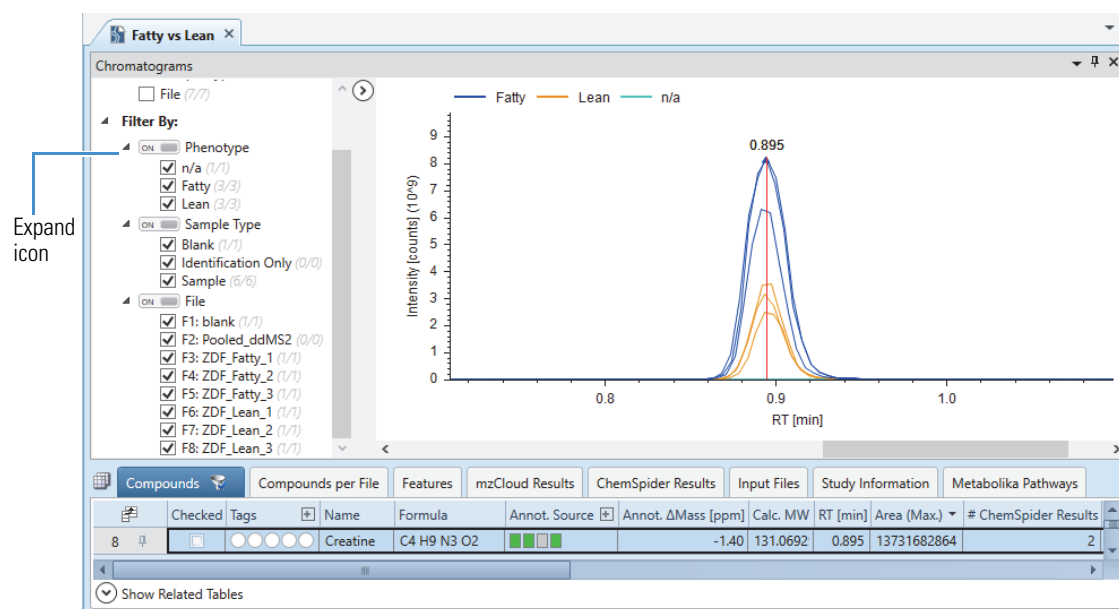
Figure 7. Overlaid traces without shading



Filter the chromatographic traces

To review the traces for creatine by phenotype, sample type, or individual files, do the following:

1. Under Filter By, expand Phenotype, Sample Type, and File by clicking their expand icons, ▲.



Note Under Group By, the count for Phenotype is 3 out of 3 (3/3). Under Filter By, Phenotype lists three groups, and all three groups (Lean, Fatty, and N/A) are selected. The N/A group consists of the Blank sample.

Under Filter By > File, the count for F2: Pooled_ddMS2 is (0/0). The application does not report compounds for the Identification Only sample type. It uses the fragmentation scans in this sample type for the mzCloud and mzVault searches and the Predict Composition node's Fragments matching algorithm.

Under Filter By > File, the count for the Blank is (1/1). When the processing workflow includes the Fill Gaps node, the application detects the compound (calculates a nonzero area value) in all the input files if it detects the compound in any of the input files.

2. Do any of the following under Filter By:

- To view only the overlaid traces for the Lean phenotype, under Phenotype, click the **N/A** and **Fatty** check boxes.
- To view only the overlaid traces for the Fatty phenotype, clear the **N/A** and **Lean** check boxes.

Review the full scan MS1 spectra

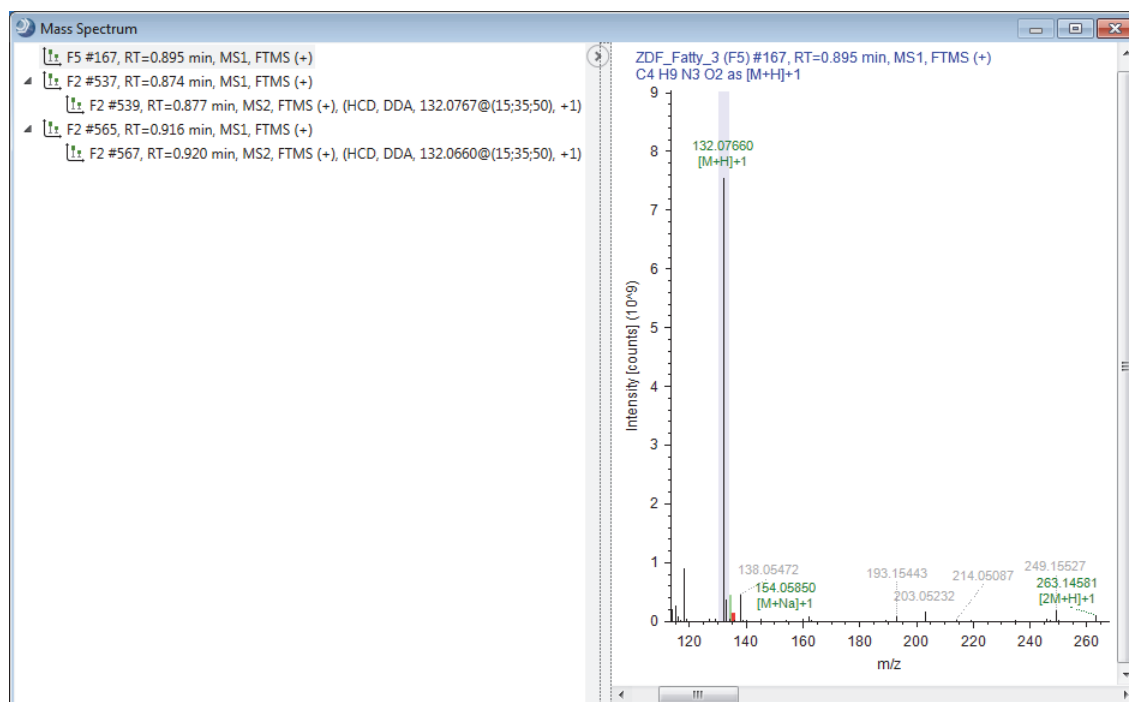
- To view only the trace for the blank sample, under Sample Type, clear the **Identification** and **Sample** check boxes. Or, clear the Lean and Fatty check boxes under Phenotype.
- To view only the trace for a specific file, under File, clear all the check boxes except for the check box for the file you want to view.

When the application predicts the formula for a compound, it annotates the isotopic peaks in the full MS1 scans.





To view the color-coded isotopes for a compound with a predicted formula in the Mass Spectrum view, do the following:

- With the Compounds table sorted in descending order by the Area (Max.) column, select **row 8** (creatine, C₄H₉N₃O₂).
- Expand the Mass Spectrum view to the full screen width by closing the Chromatograms view or by dragging the Mass Spectrum view by its title bar to a second monitor.

To close the Chromatograms view, click its **Close** icon.



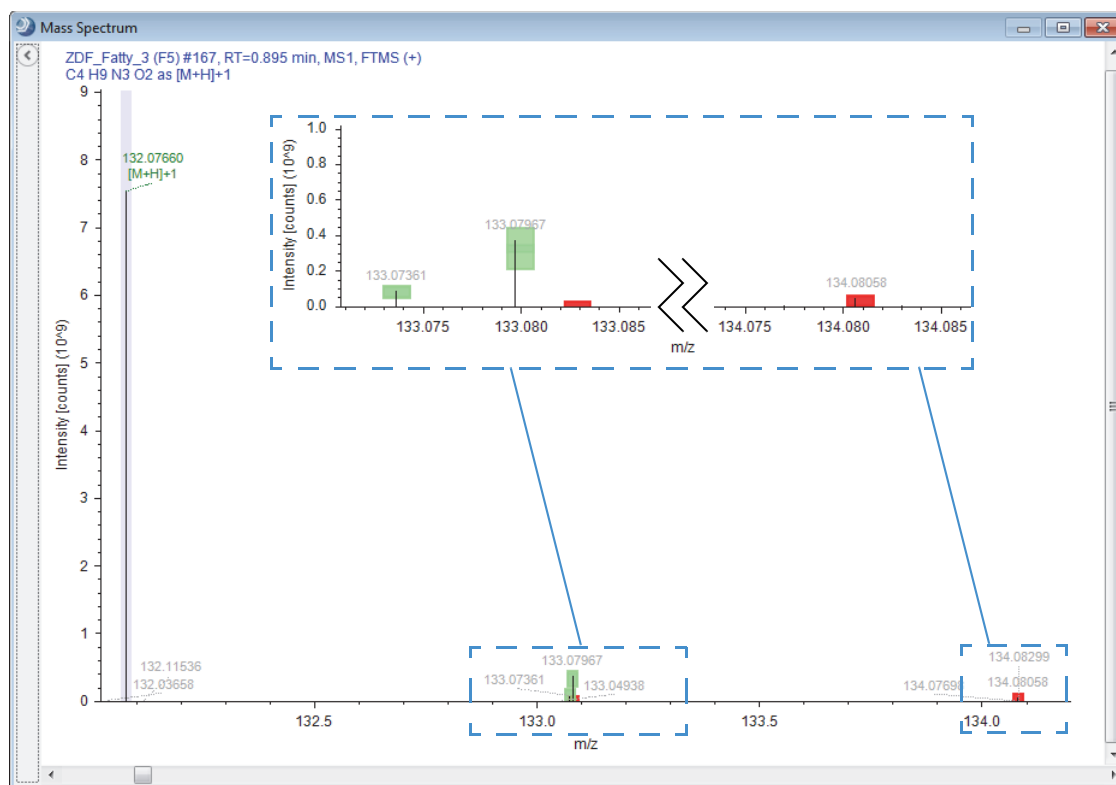
The mass spectral peaks in the full MS scan are color-coded as follows:

-  Lavender bars indicate centroids for A0 (monoisotopic) ions. The x-axis position and the width of the bar reflect the expected m/z value of the centroid and the user-specified mass tolerance, respectively.
-  Green rectangles indicate matching centroids for isotopic ions. When you zoom in on the matching centroid, the x-axis position and width of the rectangle reflect the expected m/z value of the centroid and the user-specified mass tolerance, respectively. The y-axis position and height of the rectangle reflect the expected relative intensity of the centroid and the user-specified intensity tolerance, respectively.
-  Red rectangles indicate centroids that are missing from the expected isotopic pattern.
-  Blue rectangles indicate centroids that are missing from the expected isotopic pattern but that are also expected to have an intensity below the measured baseline noise (determined by the Fourier transform mass spectrometry (FTMS) mass analyzer).

- To zoom in on the isotopes detected for the $[M+H]^+1$ adduct ion, drag the mouse pointer across the x axis from m/z 132 to m/z 135.

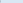


This figure shows three matching isotopes and two missing isotopes for the $[M+H]^+1$ adduct ion of creatine.

Matching	Missing
Monoisotopic A0 ion (C ₄ H ₉ N ₃ O ₂) at m/z 132.07660	A1 ion (C ₄ [2]H H ₈ N ₃ O ₂) at m/z 133.08281
A1 ion (C ₄ H ₉ [15]N N ₂ O ₂) at m/z 133.07361	A2 ([13]C ₂ C ₂ H ₉ N ₃ O ₂) ion at m/z 134.08085
A1 ion ([13]C C ₃ H ₉ N ₃ O ₂) at m/z 133.07967	



- For more information about the matching and missing isotopes, click **Show Related Tables** below the Compounds table. Then, click the **Predicted Compositions** tab.

This figure shows that the Predict Compositions node matched 3 out of 5 isotopes to the theoretical isotope pattern for the calculated formula.

Hide Related Tables														
Structure Proposals		Compounds per File		Predicted Compositions	Metabolika Results		mzCloud Results		ChemSpider Results		Metabolika Pathways			
	Checked	Compound Match	Formula	Molecular Weight	Δ Mass [Da]	Δ Mass [ppm]	RDBE	H/C	Rank	# Matched Iso.	# Missed Iso.	# Matched Frag.	SFit [%]	
1			C4 H9 N3 O2	131.06948	-0.00018	-1.40	2.0	2.2	1	3	2	8	34	

Review the fragmentation spectra

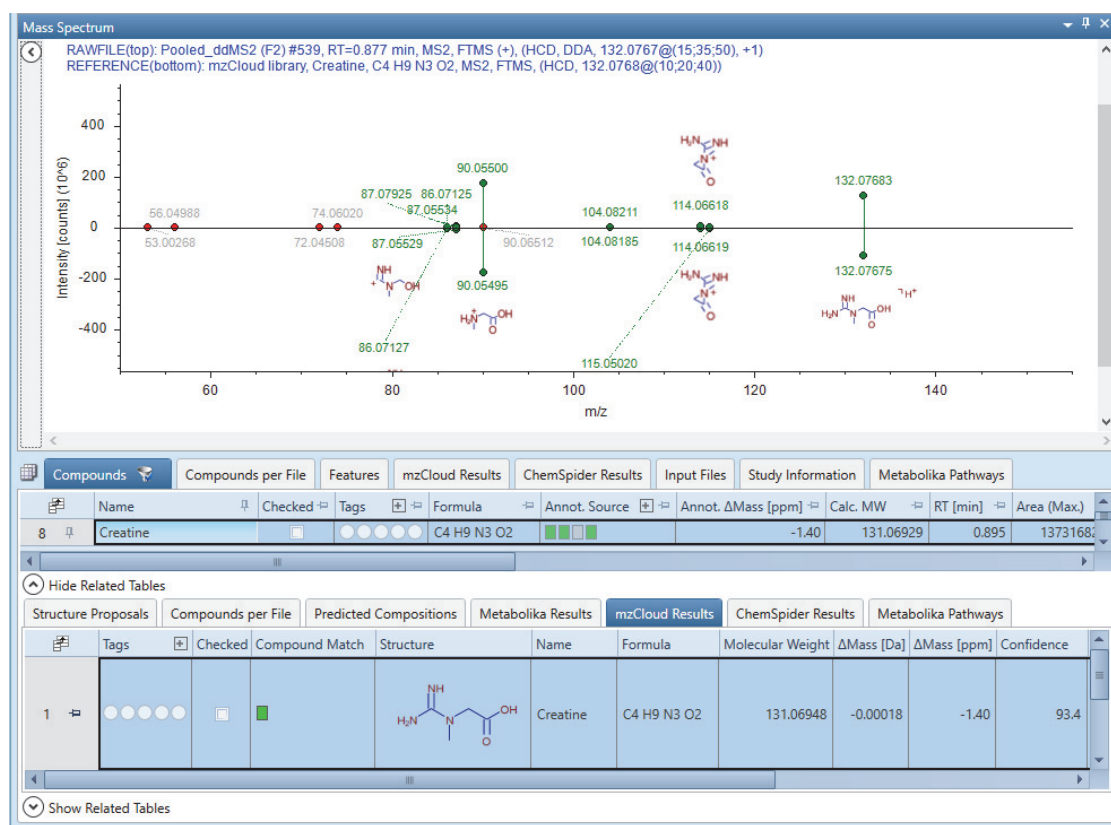
To view the results of an mzCloud search for a compound and display a mirror plot, do the following:

- With the Compounds table sorted in descending order by the Area (Max.) column, select **row 8** (Creatine, C₄ H₉ N₃ O₂).
- Click **Show Related Tables** to display the related tables below the main tables.
- Click the **mzCloud Results** tab to display the mzCloud Results table.
- Select the **row 1** in the mzCloud Results related table.

An annotated mirror plot appears in the Mass Spectrum view with the mass spectrum from the Pooled_ddMS2 sample on the top and the mzCloud reference spectrum on the bottom (Figure 8 on page 25).

A red circle on the m/z axis indicates an unmatched fragment. The green circles at the ends of centroids indicate matched fragments.

Figure 8. Annotated mirror plot for creatine



Apply the Statistics layout

The application comes with the factory default layout and four named layouts: Identification, Quantification, Stable Isotope Labeling, and Statistics. When running statistical analyses, for ease of use, apply the Statistics layout.

To apply the Statistics layout, do the following:

1. If the result file is not the active page in the application window, click the result file tab.
2. From the menu bar, choose **Window > Reset Layout**. Then, choose **Window > Apply Layout > Statistics**.

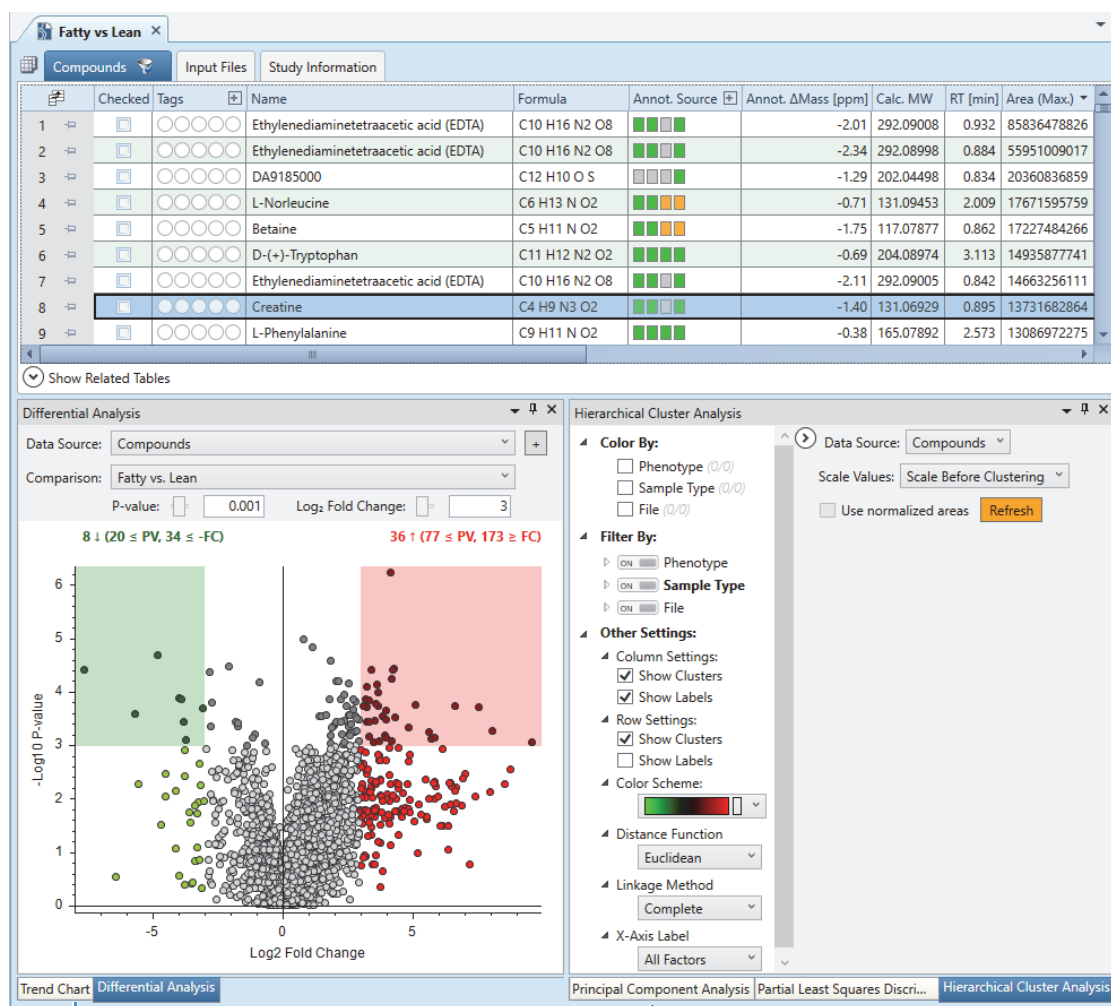
For the example result file, the Statistics layout does the following (see Figure 9):

- Closes the Chromatogram view and the Mass Spectrum view. Moves the result tables to the top of the page and hides most of the main tables, except for the Compounds table and Input Files table.
- Opens the Differential Analysis and Trend Chart views as tabbed views on the bottom left of the page with the Differential Analysis view selected as the active tab. For the differential analysis, sets the P-value parameter to 0.001 and the Log₂ Fold Change parameter to 3.
- Opens the Principal Component Analysis, Partial Least Squares Discriminant Analysis, and Hierarchical Cluster Analysis views as tabbed views on the bottom right of the page.

Note The Compounds table, the Partial Least Squares Discriminant Analysis view, the Differential Analysis view, the Hierarchical Cluster Analysis view, and the Loadings Plot page of the Principal Component Analysis view are interactive—that is, checking the points in any of these views checks the points in all of these views.

In the Hierarchical Cluster Analysis view, a segmented bar—red for checked and gray for unchecked—to the right of the heat map indicates whether a compound is checked. In the other views, checked data points turn blue.

Figure 9. Statistics layout for the example result file



Left pane

Right pane

Go to the next topic to [“View a trend chart for a single compound or a set of trend lines for multiple compounds.”](#)

View a trend chart for a single compound or a set of trend lines for multiple compounds

Use the Trend Chart view to compare the distribution of the chromatographic peak areas for each compound by sample group. When you select a single compound in the Compounds table, you can view the distribution as a box-and-whiskers plot or as a trend line plot. When you select multiple compounds in the Compounds table, the application automatically displays the distribution for each compound as a trend line plot.

To learn how to work with the Trend Chart view, follow these topics:

- [View a box-and-whiskers plot for a single compound](#)
- [View the trend line plots for multiple compounds](#)

View a box-and-whiskers plot for a single compound

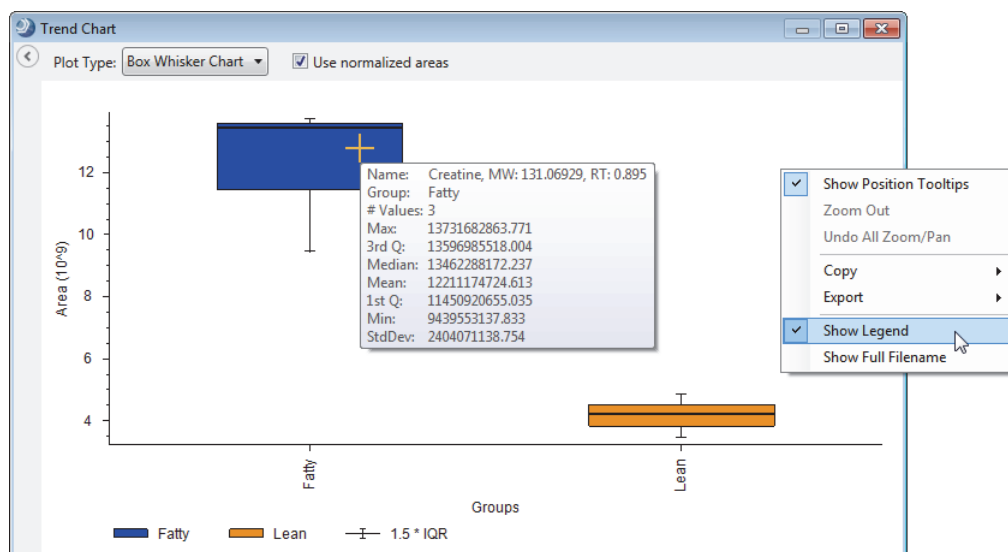
To view a box-and-whiskers plot for a single compound, do the following:

1. Apply the Statistics layout (see [“Apply the Statistics layout”](#) on page 25).

Note Applying the Statistics Layout changes the sort order of the Compounds table.

2. Sort the Compounds table in descending order by Area (Max.).
3. Click the **Trend Chart** tab in the left pane.
4. Select **row 8** (Creatine, C4 H9 N3 O2).
5. Right-click the chart and choose **Show Legend**.

6. Place the cross-hair cursor anywhere on a box or whisker to display a tooltip with descriptive statistics. This figure shows the box-and-whiskers plot for creatine and the Tooltip for the Fatty sample group.

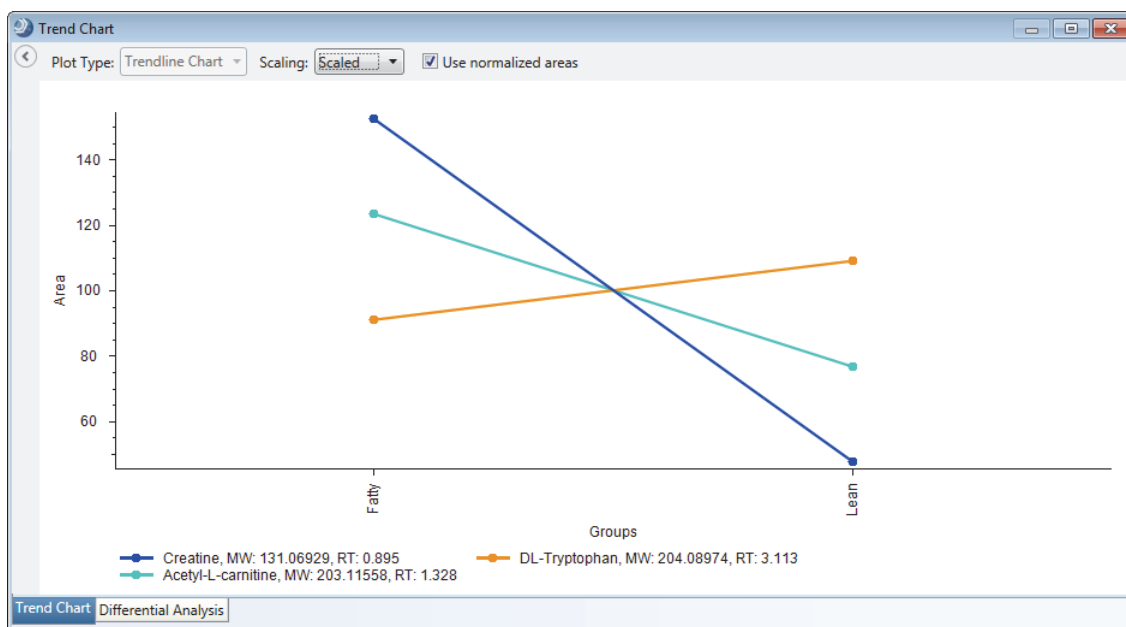


View the trend line plots for multiple compounds

To view the trend line plots for multiple compounds, do the following:

1. Hold down the CTRL key and select the following rows in the Compounds table: **row 8** (creatine), **row 6** (D-(+)-tryptophan), and **row 14** (acetyl-L-carnitine).

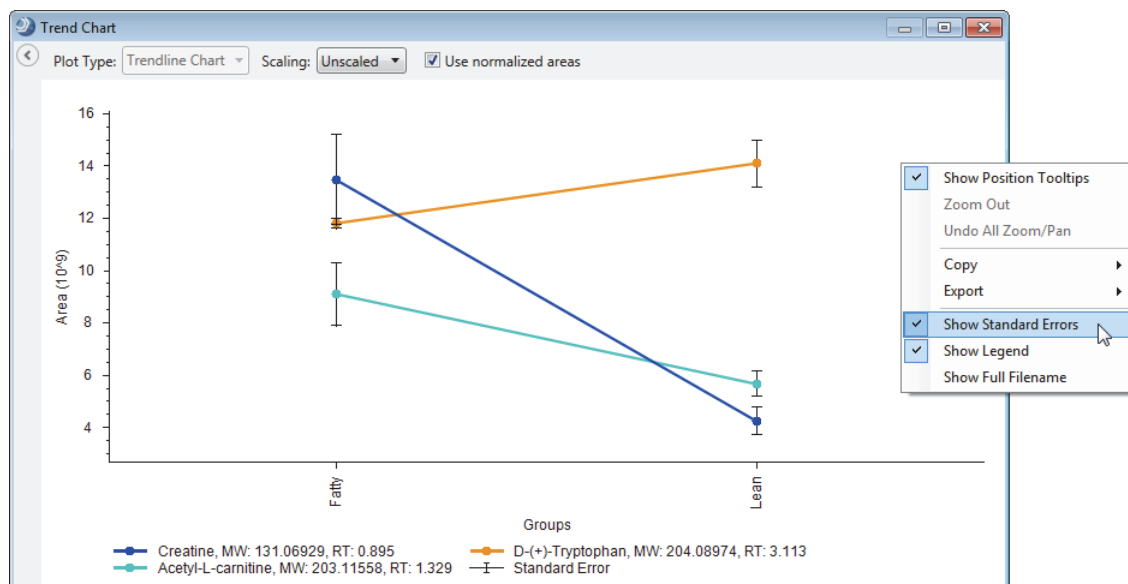
This figure shows the trendline plots for the selected compounds in the Trend Chart view.



2. To display the error bars for the trend line plots, do the following:
 - a. In the Scaling list, select **Unscaled**.
 - b. Right-click the plot and choose **Show Standard Errors**.

Visualize the detected compounds as a volcano plot

This figure shows trend line plots with error bars for the standard error.



Use the Differential Analysis view to select the significant compounds for the comparison groups, which in this case are the Fatty and Lean ZDF rats.

The volcano plot is a plot of the p-value, the result of a significance test, on the y axis versus the log₂ fold change between two sample groups on the x axis. The y-axis scale is the $-\log_{10}$ of the p-value. As the p-value increases from 0 to 1, the $-\log_{10}$ of the p-value decreases from infinity to 0 ($-\log_{10} 0.05$ equals 1.3).

The compounds in the shaded regions of the volcano plot are statistically significant for the specified p-value and outside the specified upper and lower fold change thresholds. The compounds in the green-shaded area are down-regulated compounds, and the compounds in the pink-shaded area are up-regulated compounds.

If you have not already applied the Statistics Layout, apply it now (see [“Apply the Statistics layout”](#) on page 25). Then, sort the Compounds table by the Area (Max) column.

To view the result of the differential analysis that you set up on the Groups and Ratios page, do the following:

1. Click the **Differential Analysis** tab.
2. To display the legend for the volcano plot, right-click the view and choose **Show Legend**.

The legend appears at the bottom of the page. Each data point represents a compound. Placing the crosshair cursor over a data point displays a tooltip with the XY coordinates, MW, RT, maximum area, and number of adducts.

Figure 10 shows the default Statistics layout for the Differential Analysis view, the added legend, the tooltip for the data points, and the shortcut menu for the view

Figure 10. Default settings for the Statistics layout



3. To learn how to work interactively with the Compounds table and the Differential Analysis view, do the following:
 - a. To select the up-regulated compounds or the down-regulated compounds, right-click the plot and choose **Check All Up-regulated Points**.
The checked data points turn blue, and the application checks the corresponding check boxes for the selected data points in the Compounds table.



- b. Right-click the Differential Analysis view and choose **Uncheck All Visible Data Points**.

c. To view information about the compound for a data point in the Compounds table, do one of the following:

- Double-click the data point in the Differential Analysis view.

The application automatically scrolls to the selected compound in the Compounds table, but it does not select the check box for the compound.

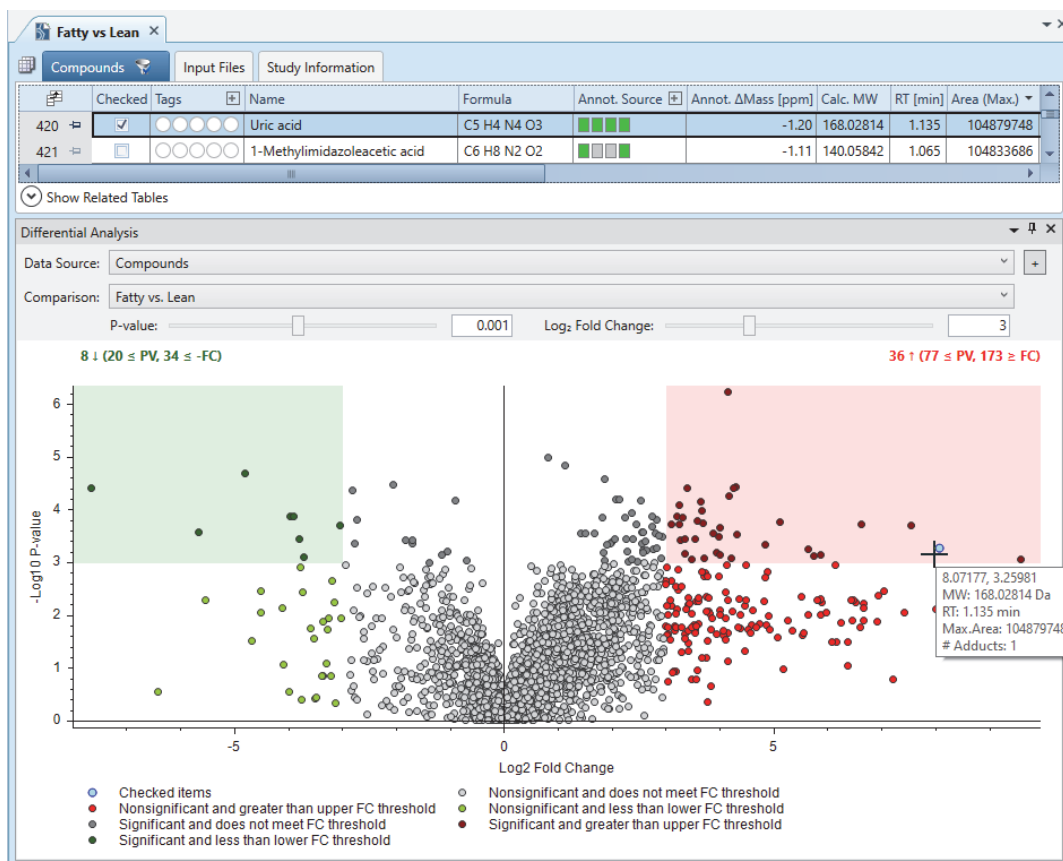
—or—

- Right-click the data point and choose **Check Point**.

The data point turns blue (Figure 11). The application automatically scrolls to the selected compound in the Compounds table and selects the check box for the compound.

d. Right-click the Differential Analysis view and choose **Uncheck Point**.

Figure 11. Checked data point



View the principal component analysis

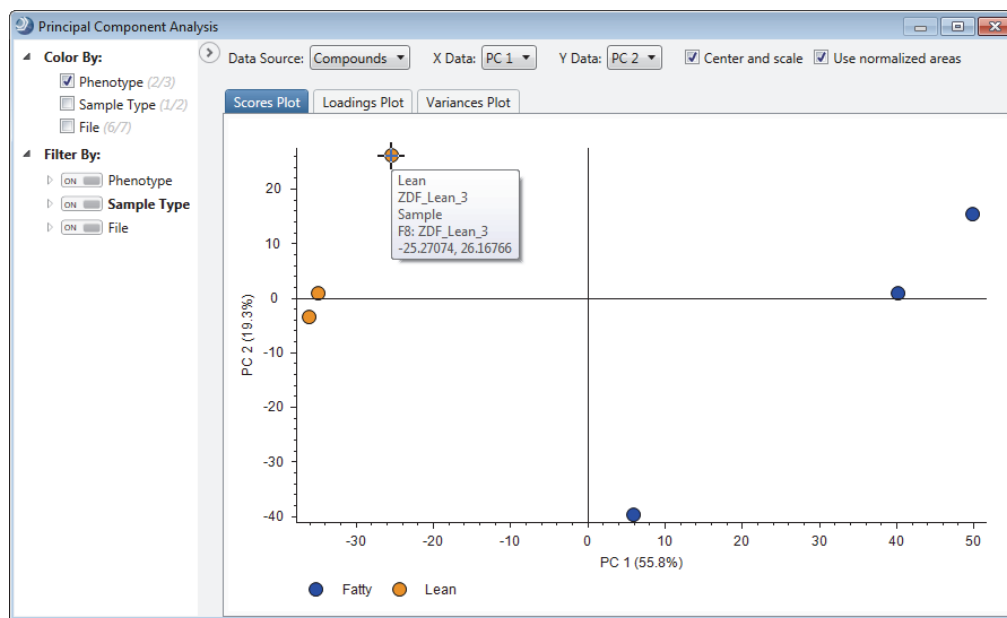
Use the Principal Component Analysis view to display the results of the principal component analysis. The principal component analysis reduces the dimensionality of the data set to a set of principal components, PC1, PC2, and so on, where PC1 is the principal component with the most variance.

To view the principal component analysis, do the following:

1. Apply the Statistics layout (see [“Apply the Statistics layout”](#) on page 25).
2. Click the **Principal Component Analysis** tab in the right pane.

Use the Scores Plot page to interpret the relationship among the sample groups. Sample groups that are near each other are similar.

This figure shows the Scores Plot page for the Fatty and Lean sample groups.



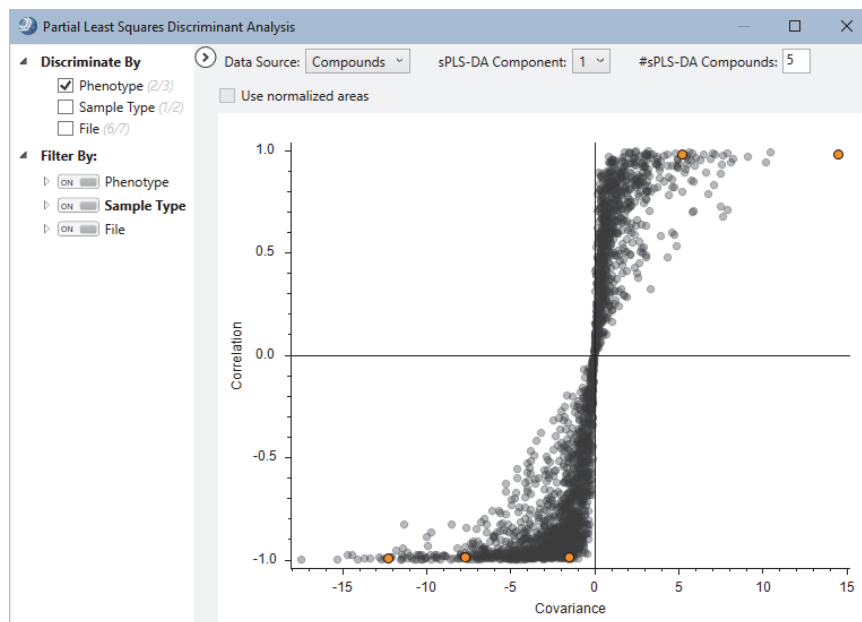
Find a set of discriminating compounds

Use the Partial Least Squares Discriminant Analysis view to find a set of compounds that you can use to discriminate between the two sample groups.

To find a set of discriminating compounds, do the following:

1. Apply the Statistics layout (see [“Apply the Statistics layout”](#) on [page 25](#)).
2. Click the **Partial Least Squares Discriminant Analysis** tab in the right pane.
3. If the graph includes checked data points (highlighted in blue), right-click the view and choose **Uncheck All Visible Points**.

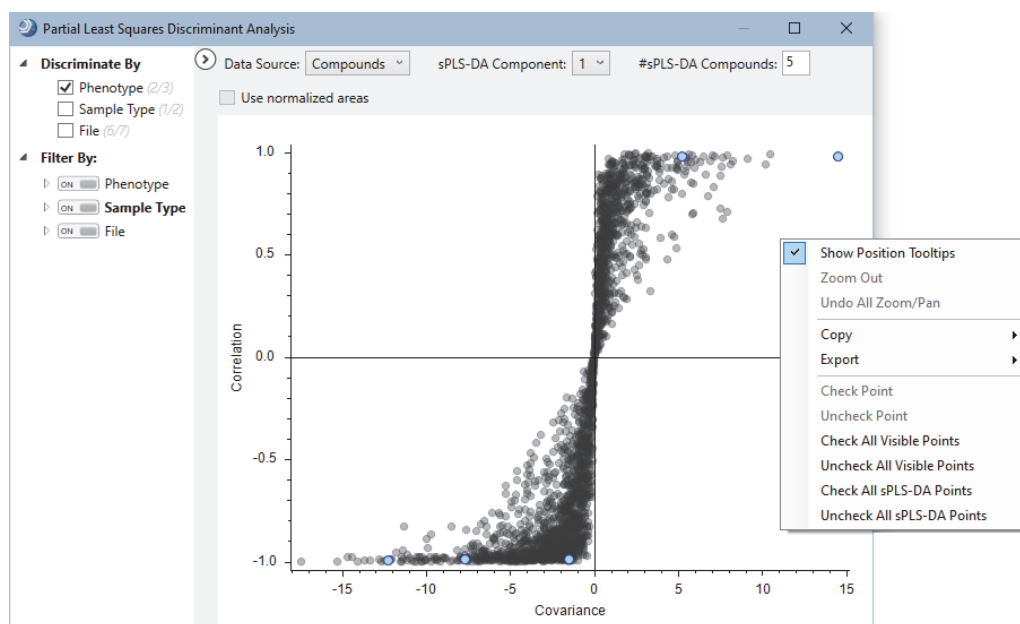
This figure shows the Partial Least Squares Discriminant Analysis view with the default settings for the Statistics layout. The orange circles represent the five discriminating compounds for the two phenotypes.



- To automatically select the check boxes for the discriminating compounds in the Compounds table, right-click the view and choose **Check All sPLS-DA Points**.

The color of the circles changes to blue and the check boxes for the associated compounds in the Compounds table are selected.

Figure 12. Five selected compounds



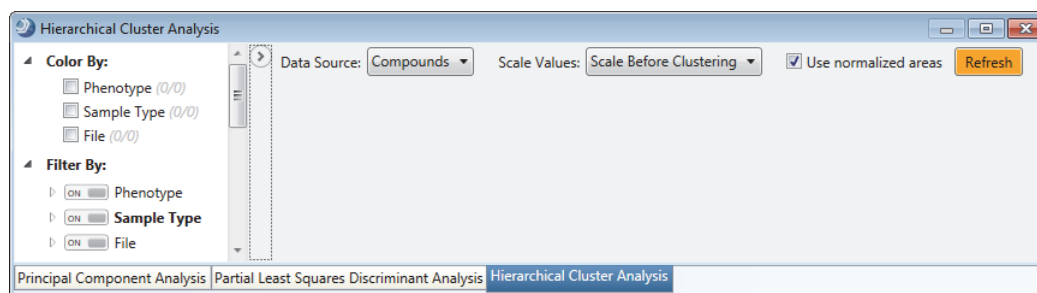
Tip If the discriminating compounds are checked, right-click the view and choose **Uncheck All sPLS-DA Points** before changing the number of discriminating compounds.

To display only the checked compounds in the Compounds table, see “[Filter the compounds their checked status](#)” on [page 38](#).

Use the Hierarchical Cluster Analysis view to visualize the correlation between detected compounds and selected samples in a two-dimensional array of color-coded rectangles (heat map) where each rectangle represents the relative amount (by area) of a specific compound in a specific sample.

To run a hierarchical cluster analysis for the compounds in the Compounds table, do the following:

- Apply the Statistics layout (see “[Apply the Statistics layout](#)” on [page 25](#)).
- Click the **Hierarchical Cluster Analysis** tab in the right pane.



- In the Hierarchical Cluster Analysis view, click **Refresh**.

A heat map and two dendograms for the cluster analysis appear.

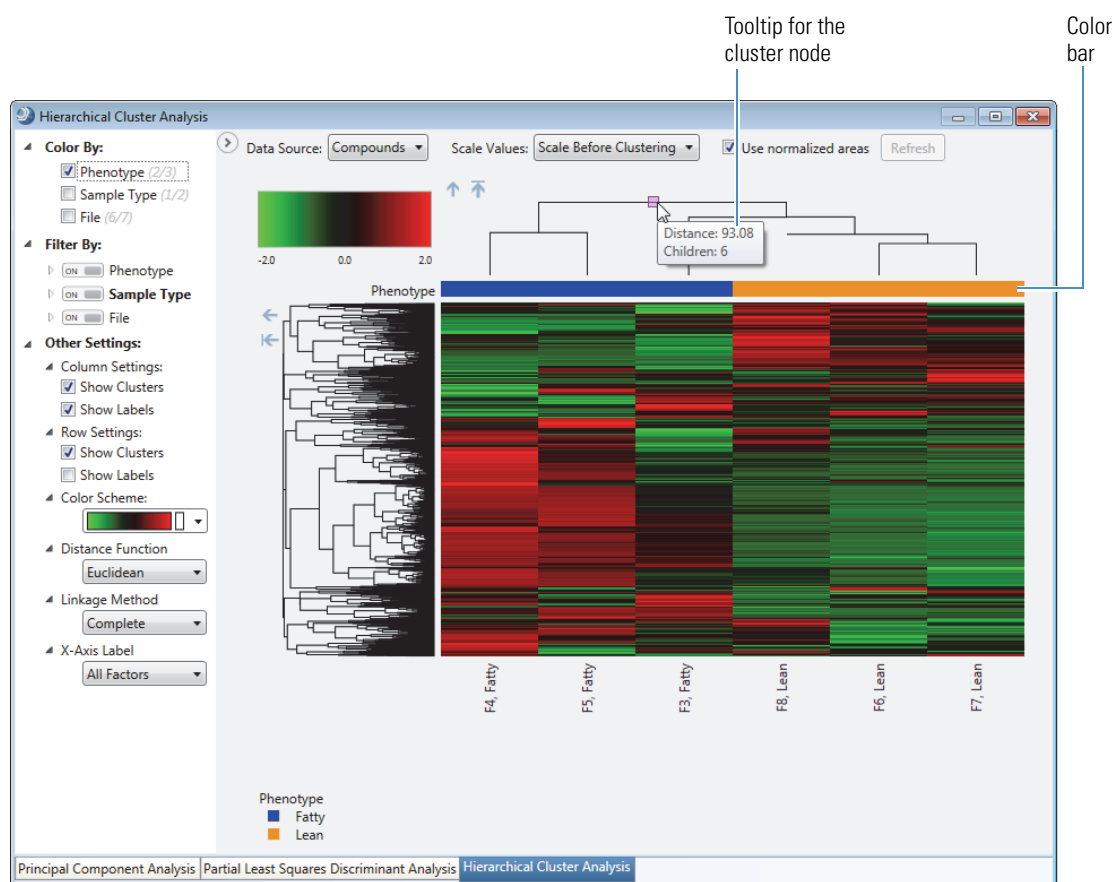
- In the left pane, under Color By, select the **Phenotype** check box.

A color bar appears above the heat map to visually differentiate the samples by the two phenotypes—Fatty and Lean ([Figure 13](#)).

Run a hierarchical clustering analysis

Figure 13 shows a hierarchical cluster analysis for the compounds in the Compounds table. Pointing to a cluster's node displays a tooltip.

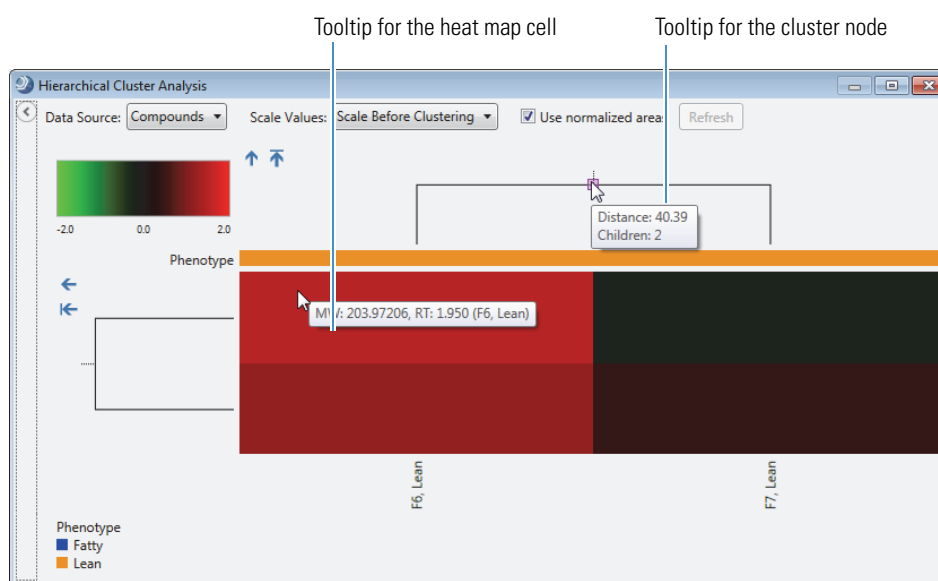
Figure 13. Heat map



5. To zoom in on a specific area of the heat map, drag the pointer across the rectangular area of interest.
6. To view information for a cell in the heat map, point to it.

The tooltip displays the row and column coordinates. The row coordinates are the compound's name, MW, and RT. The column coordinates are the file ID and study factor values.

This figure shows a magnified view of four cells in the upper right corner of the heat map and their corresponding dendograms.



View the Metabolika pathways for a compound

The Map to Metabolika Pathways node searches for compounds that match the detected compounds in the selected Metabolika Pathway files.

To view a Metabolika pathway for a detected compound, do the following:

1. From the menu bar, choose **Window > Reset Layout**. Then, at the prompt, click **Yes**.
2. Sort the Compounds table by the Area (Max.) column in descending order.
3. Select **row 10** (L-Tyrosine).
4. Below the Compounds table, click **Show Related Tables**.
5. Click the **Metabolika Pathways** tab.
6. In the Metabolika Pathways table, select **row 5** (L-tyrosine degradation IV (to 4-methylphenol)).

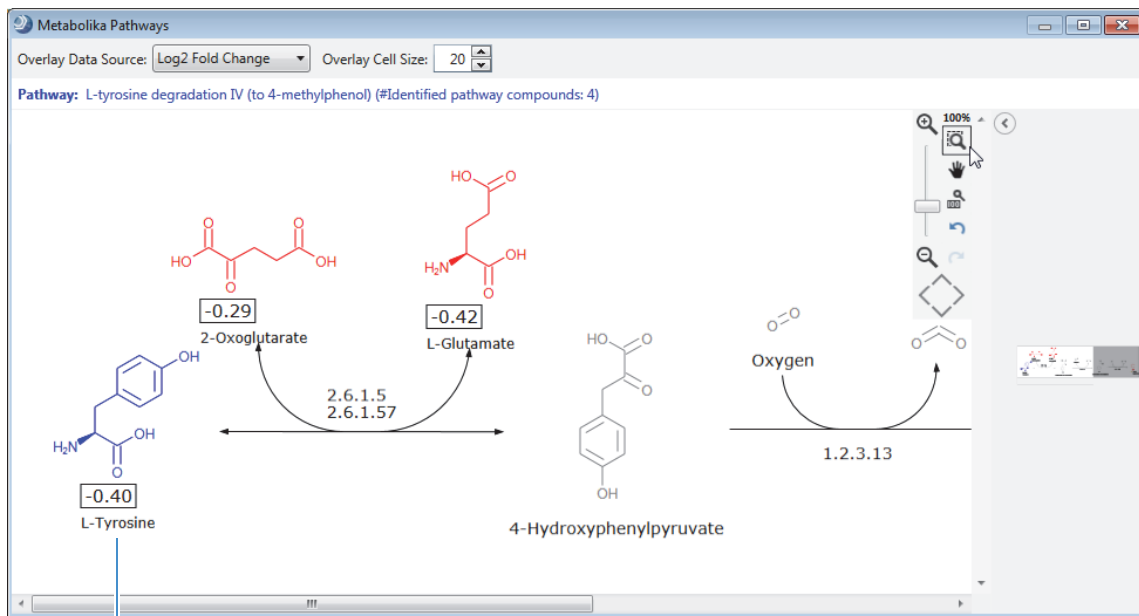
#	Checked	Tags	Pathway Name	Metabolika Compound Ids	Metabolika Compound Names	Metabolika Compound Formula	# Mapped Compounds	# Matched Compounds	# Compounds in Pathway
1	<input type="checkbox"/>		Superpathway of chorismate metabolism	692	L-Tyrosine	C9 H11 N O3	17	15	152
2	<input type="checkbox"/>		Superpathway of aromatic amino acid biosynthesis	692	L-Tyrosine	C9 H11 N O3	15	12	47
3	<input type="checkbox"/>		Superpathway of thiamine diphosphate biosynthesis I	692	L-Tyrosine	C9 H11 N O3	11	6	44
4	<input type="checkbox"/>		Superpathway of betaine biosynthesis	692	L-Tyrosine	C9 H11 N O3	10	8	59
5	<input checked="" type="checkbox"/>		L-tyrosine degradation IV (to 4-methylphenol)	692	L-Tyrosine	C9 H11 N O3	9	6	9
6	<input type="checkbox"/>		Superpathway of 4-hydroxybenzoate biosynthesis (yeast)	692	L-Tyrosine	C9 H11 N O3	8	7	22
7	<input type="checkbox"/>		Superpathway of L-tyrosine biosynthesis	692	L-Tyrosine	C9 H11 N O3	7	5	28
8	<input type="checkbox"/>		Superpathway of rosmarinic acid biosynthesis	692	L-Tyrosine	C9 H11 N O3	7	6	37
9	<input type="checkbox"/>		Novobiocin biosynthesis	692	L-Tyrosine	C9 H11 N O3	6	5	60
10	<input type="checkbox"/>		Superpathway of plastoquinol biosynthesis	692	L-Tyrosine	C9 H11 N O3	6	5	17
11	<input type="checkbox"/>		Superpathway of L-phenylalanine and L-tyrosine biosynthesis	692	L-Tyrosine	C9 H11 N O3	5	4	11

7. From the application menu bar, choose **View > Metabolika Pathways**.

The view opens to the right of the result tables and displays the selected pathway. The structure for the selected compound is blue, the structures for other compounds in the Compounds table are red, and the structures for other compounds in the pathway that are not in the Compounds table are black.

8. In the Overlay Data Source list, select **Log2 Fold Change**.
9. In the Overlay Cell Size list, increase the value to **20**.

This figure shows the selected pathway, data source, and overlay cell size.



Compound selected in the Compounds table

Review the results of the mzLogic analysis

An mzLogic analysis requires candidate structures from the ChemSpider database, a Metabolika pathway, or a mass list with structures. It also requires matching structures from an mzCloud similarity search.

Tip If the processing workflow for an analysis did not include the Search ChemSpider node or the Apply mzLogic node, you can run an mzLogic Analysis from the mzLogic Analysis view.

To review the results of an mzLogic analysis for structure candidates, do the following:

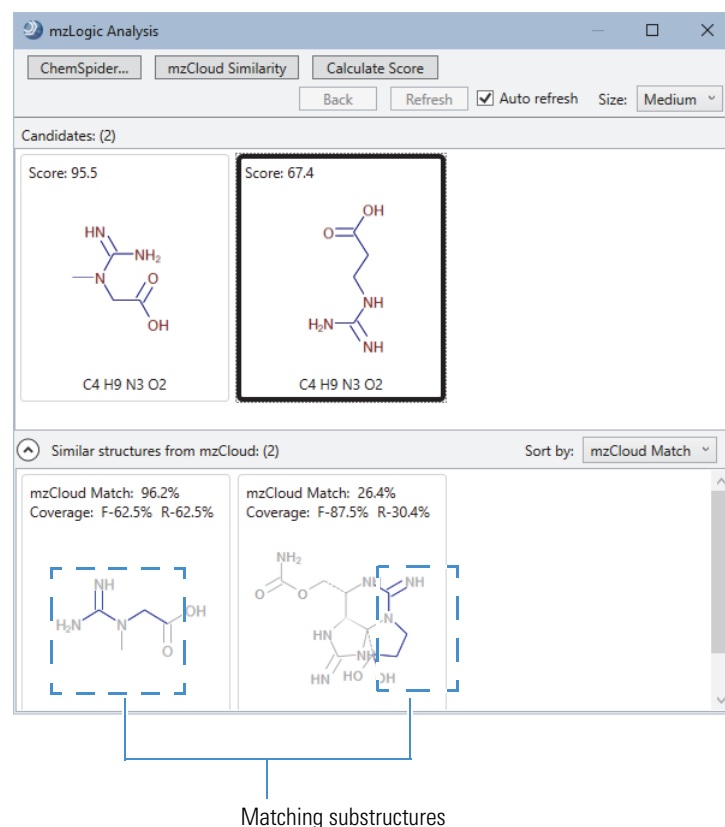
1. In the Compounds table (sorted in descending order by the Area (Max.) column, select **row 8** (creatine).
2. From the application menu bar, choose **View > mzLogic Analysis**.

The mzLogic Analysis view opens to the right of the result tables.

3. Below the Candidates pane, click **Similar Structures from mzCloud**.
4. In the Candidates pane, select the structure for the lower ranking hit with a score of 67.4.

Figure 14 shows the matching substructure in blue for the lower ranking compound in the related ChemSpider Results table for creatine.

Figure 14. mzLogic Analysis results



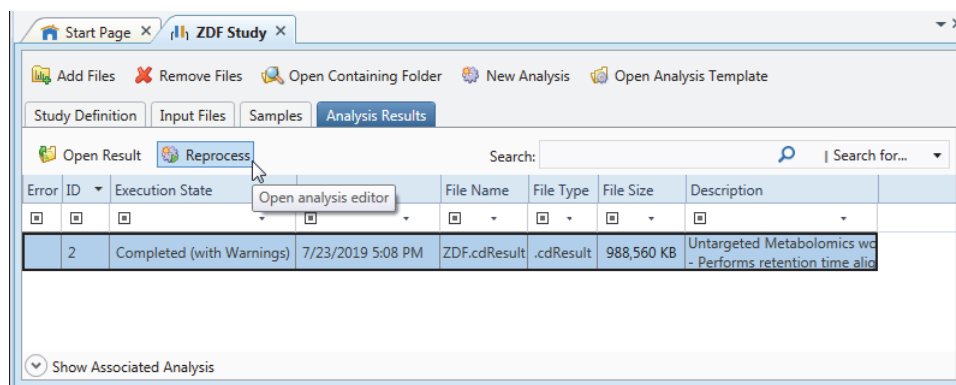
Modify the processing workflow and partially reprocess the analysis

You can modify the parameter settings for any of the scoring, mapping, or search nodes or add or remove these nodes from the processing workflow without reprocessing the entire workflow—that is, without reprocessing the core nodes.

Note In the processing workflow for this tutorial, the Input Files, Select Spectra, Align Retention Times, Detect Compounds, Group Compounds, and Fill Gaps nodes are core nodes.

To partially reprocess the same set of input files with a modified processing workflow, do the following:

1. Open the Analysis Results page of the study.
2. Select the analysis that you want to reprocess.



3. Click **Reprocess**.

The Grouping & Ratios and Workflows tabs appear, and a Caution symbol appears in the Analysis pane to the right of the Edit button. Pointing to the Caution symbol displays the following message:

Unchanged workflow. No node chosen for processing.

4. Click the **Workflows** tab.

All the nodes have lost their distinguishing colors, and the right sides of all the nodes have a gray tab.

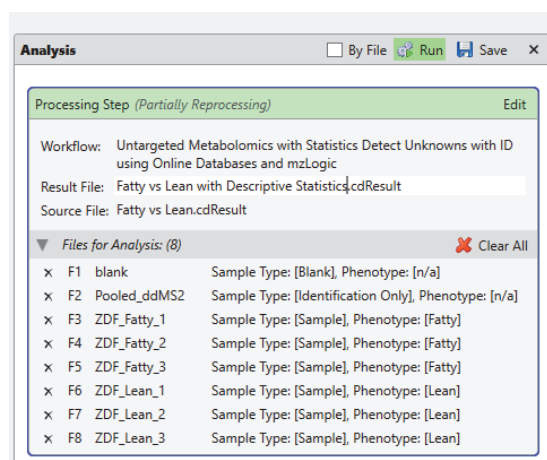
5. Add or remove any of the pathway mapping, compound scoring, or post-processing nodes.

For example, add the Descriptive Statistics node by dragging it from the Workflow Nodes pane to the Post-Processing Nodes pane at the bottom of the Workflows page.

When you modify the downstream nodes, the application creates a Source file box where it copies the file name of the original file in the Analysis pane. You cannot edit the original file name.

6. In the Result File box, rename the result file that you want the application to create by partially reprocessing the data.

Figure 15. Analysis pane set for partial reprocessing



7. Close the original result file.

8. Click **Run**.

One of these actions occurs:

- If the original result file is still open, the application prompts you to close it. At the prompt, click **Abort**, close the original result file, and then click **Run**.
- If the original result file is closed, the application submits the run to the job queue. Reprocessing the data after adding the Descriptive Statistics node takes less than a minute.

Filter the data and export the analysis results

Use the result filters to select the compounds of interest

To create a report for your records or a file for external use by the TraceFinder™ application or the Xcalibur data system, filter the compounds table to display only the compounds of interest, and then export the results using the appropriate format.

To filter the compounds table and export the results, do the following:

1. [Use the result filters to select the compounds of interest](#)
2. [Export the results to a spreadsheet, text, or CSV file](#)

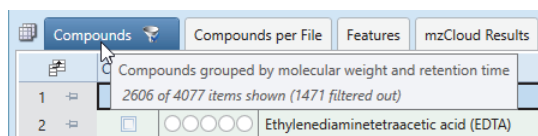
The analysis found a total of 4077 compounds, including 1471 background compounds. By default, background compounds are hidden, so the Compounds table lists 2606 compounds.

To reduce the number of compounds to review—for example, to display only compounds that are statistically significant, filter the table by p-value. To reduce the number of compounds to the discriminating compounds for the comparison groups or to the compounds of interest for other reasons, check the discriminating compounds or the compounds of interest to you and filter the table by the checked property.

For details, see these topics:

- [Filter the compounds by their p-values](#)
- [Filter the compounds their checked status](#)

Note Pointing to the Compounds tab or the scroll bar displays a tooltip with the number of compounds in the table.

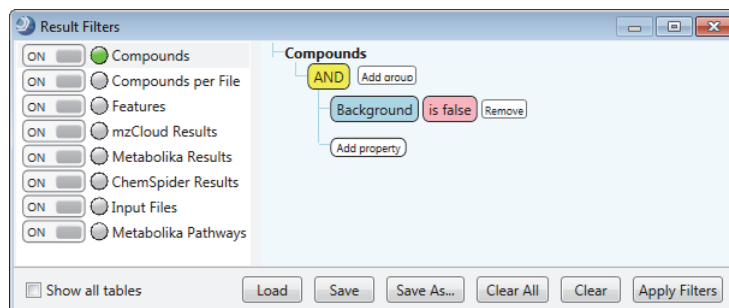


Filter the compounds by their p-values

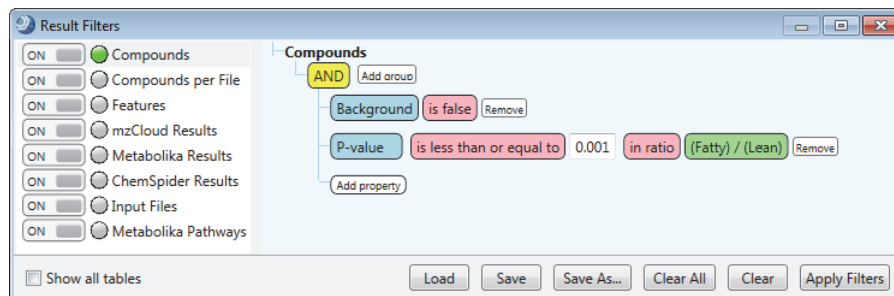
To reduce the number of compounds by using the p-value as a filter, do the following:

1. In the main tables pane, click the **Compounds** tab.
2. From the menu bar, choose **View > Result Filters**.

The Result Filters view opens as a floating window. Because the processing workflow included the Mark Background Compounds node, the filter for the Compounds table already includes a Background filter and the table does not display the background compounds.



3. In the right pane of the Result Filters view, set up a filter for the p-value as follows:
 - a. Click **Add Property** and select **P-Value** from the list.
 - b. In the pink relation list, select **Is Less Than or Equal To**.
 - c. In the value box next to the relation list, type **0.001**.
 - d. In the pink condition list, select **In Ratio**.
 - e. In the Green sample list, select **Fatty/Lean**.



4. Click **Apply Filters**.

The filter reduces the number of displayed rows in the Compounds table to 125.

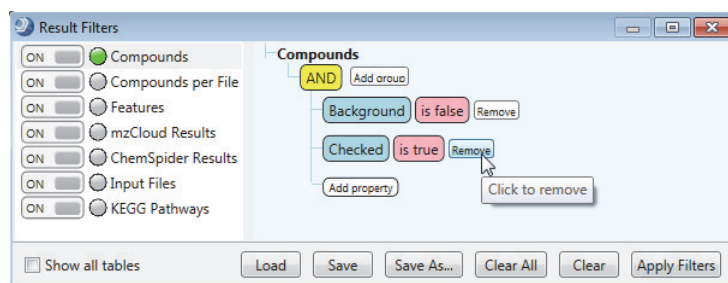
Filter the compounds their checked status

To filter the Compounds table by the Checked column, do the following:

1. If the Compounds table is not the active table, click its tab to make it active.
2. Do one of the following:
 - In the Partial Least Squares Discriminant Analysis view, select a set of discriminating compounds as described in [“Find a set of discriminating compounds”](#) on page 31.
 - In the Differential Analysis view, select the compounds of interest by checking the up-regulated compounds, the down-regulated compounds, or the visible points. See [“Visualize the detected compounds as a volcano plot”](#) on page 28.
 - Manually select the check boxes for the compounds of interest.
3. From the menu bar, choose **View > Result Filters**.
- Because the processing workflow included the Mark Background Compounds node, the Compounds table is currently filtered by one property—Background.
4. If there are more filter properties, remove them by clicking **Remove** to the right of each filter.
5. Click **Add Property** and select **Checked**.
6. Click **Apply Filters**.

The Compounds table displays only the selected compounds.

This figure shows the filter tree for the Compounds table.



7. To undo the Checked filter, click **Remove** to its right. Then, click **Apply Filters**. The Compounds table contains the original set of compounds.

Export the results to a spreadsheet, text, or CSV file

Export compounds to a spreadsheet

Before exporting the results to a spreadsheet, an Xcalibur™ inclusion/exclusion list, a TraceFinder™-compatible CSV file, or a new mzVault library, filter the results table as described in [“Use the result filters to select the compounds of interest”](#) on [page 37](#).

To learn about the export feature, do the following:

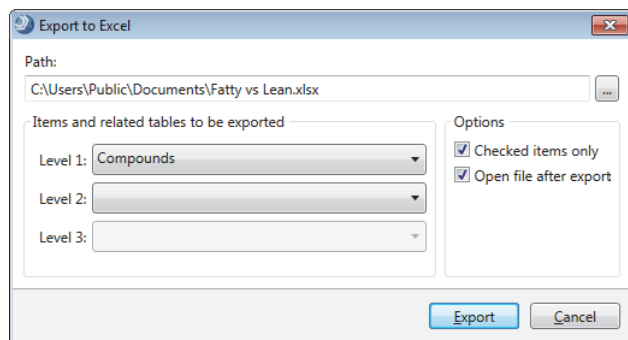
- [Export compounds to a spreadsheet](#)
- [Create an Xcalibur inclusion/exclusion list](#)
- [Create a file to import into the TraceFinder compounds database](#)
- [Export a compound to an existing mzVault library](#)
- [Export compounds to a new library](#)

To export a list of compounds to an Excel spreadsheet, do the following:

1. Do any of the following:
 - To check the number of table rows, point to the Compounds tab or the vertical scroll bar to the right of the Compounds table.
A tooltip appears with the row count.
 - To sort the table rows, click the column heading that you want to sort by once or twice to set the appropriate sort order. See [“Sort the rows in a result table”](#) on [page 20](#).
 - To display the table columns that you want to export, open the Field Chooser box and select the check boxes for the columns of interest and clear the other check boxes (see [“Show or hide table columns”](#) on [page 19](#)).
For example, to view the areas for each input file, select the **Areas** check box.
2. To export the filtered and sorted results to an Excel spreadsheet, do the following:
 - a. Right-click the Compounds table and choose **Export > As Excel**.

The Export to Excel dialog box opens.

Figure 16. Export to Excel dialog box

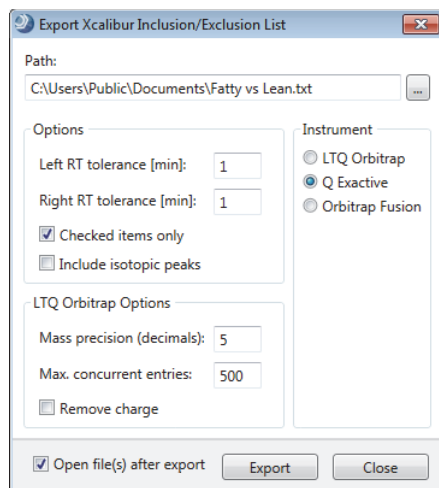


- b. Check the file name and location in the Path box. Change the file name and location as appropriate.
- c. In the Options area, select the appropriate options.
- d. Click **Export**.
- e. At the prompt, click **OK**.

Create an Xcalibur inclusion/exclusion list

To export the results to an Xcalibur inclusion/exclusion list, do the following:

1. Right-click the Compounds table and choose **Export > As Xcalibur Inclusion/Exclusion List**.

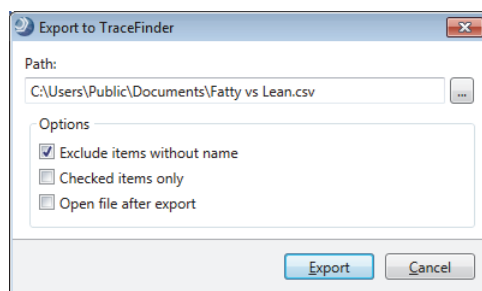


2. In the Instrument area, select the instrument format for the exported masses.
The example raw data files were acquired with a Q Exactive Orbitrap mass spectrometer.
3. (Optional) To export only the selected compounds, select the **Checked Items Only** check box.
4. (Optional) To automatically open the file in a spreadsheet application, select the **Open File(s) after Export** check box.
5. Click **Export**.
6. At the prompt, click **OK**.

Create a file to import into the TraceFinder compounds database

To export the results to a file that is compatible with the TraceFinder Compounds Database, do the following:

1. Right-click the Compounds table and choose **Export > As TraceFinder List**.
2. Select one or more of the options.

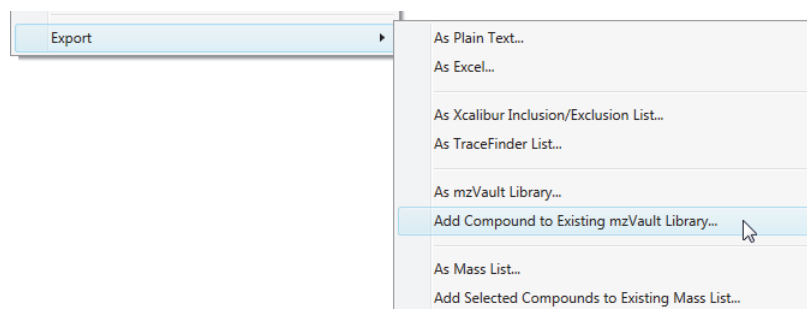


3. Click **Export**.
4. At the prompt, click **OK**.

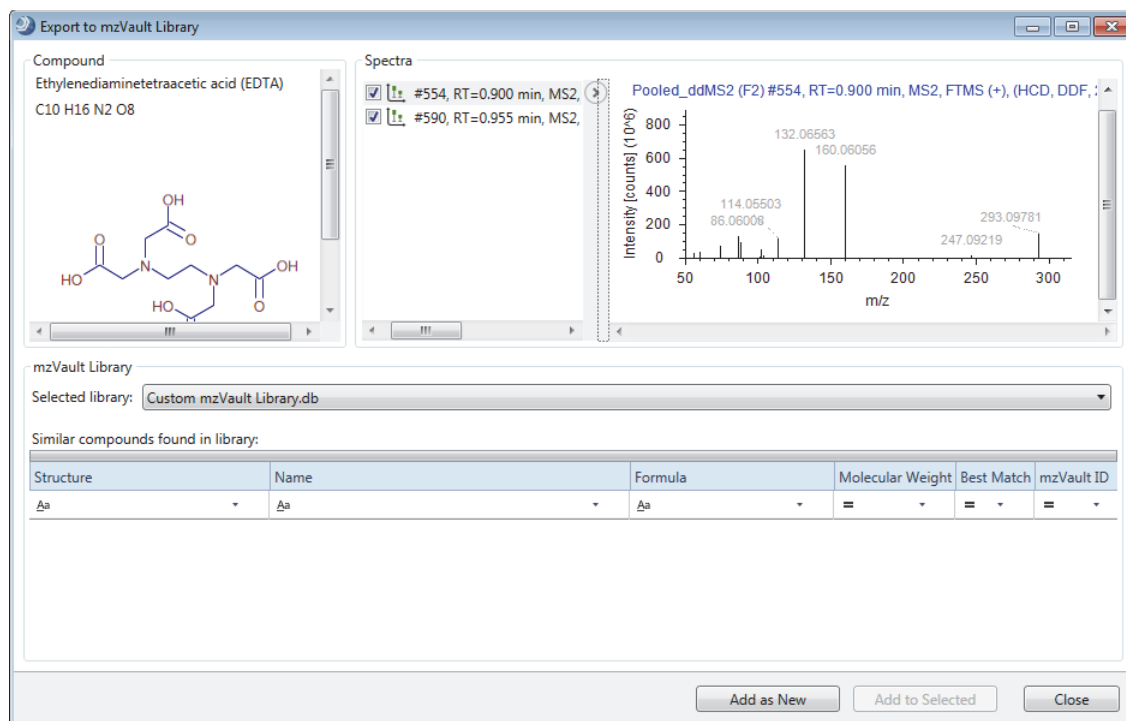
Export a compound to an existing mzVault library

To export a compound to an existing mzVault Library, do the following:

1. Right-click the compound in the Compounds table and choose **Export > Add Compound to Existing mzVault Library**.



The Export to mzVault Library dialog box opens. Its Spectra view displays the available fragmentation spectra for the selected compound.



2. Do the following:
 - In the spectrum tree, select the check boxes for the spectra that you want to add to the compound entry.
 - In the mzVault Library area, select the existing mzVault library from the Selected Library list.
3. Do one of the following:
 - To add the spectra to an existing compound entry in the library, select the entry in the Similar Compounds Found in Library list. Then, click **Add to Selected**.
 - To add the compound and the selected spectra as a new compound entry, click **Add as New**.
4. Click **Close**.

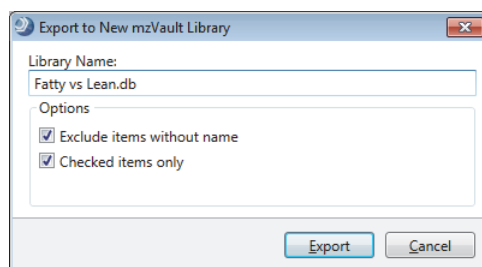
Export compounds to a new library

To export the checked compounds to a new mzVault Library, do the following:

1. Right-click the Compounds table and choose **Export > As mzVault Library**.

The Export to New mzVault Library dialog box opens. By default, the Exclude Items Without Name check box is selected.

2. Select the **Checked Items Only** check box.



3. Click **Export**.
4. At the prompt, click **OK**.

Trademarks

Compound Discoverer, Exactive Plus, FreeStyle, mzCloud, mzVault, Q Exactive, and TraceFinder are trademarks, and Exactive, Orbitrap, Orbitrap Fusion, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

Microsoft and Excel are registered trademarks of Microsoft Corporation in the United States and other countries.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.