



Agilent MassHunter Workstation Software

Report Builder Familiarization Guide



Notices

© Agilent Technologies, Inc. 2021

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

G3336-90061

Edition

November 2021

Printed in USA

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

Warranty

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

Technology Licenses

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

Restricted Rights Legend

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

Safety Notices

CAUTION

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

WARNING

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

Contents

1 Introduction

- Access the Report Builder **5**
- Access Report Builder PDF Templates **6**
- Before You Begin **7**

2 Report Template Workflows

- Quantitative Analysis - Calibration Report Template **9**
- Unknowns Analysis - Area Percent Report Template **50**

3 Modify a Report Template

- Replace the Agilent Logo with Your Own **98**
- Replace a Column in the Report **99**
- Replace a column using the Textbox Properties window **101**
- Remove a Column From the Report **103**
- Adjust the Width of a Column **104**
- Filter Results Using Special Conditions **105**
- Adjust the Size of a Graphic **107**
- Add a New Column to an Existing Report **108**
- Edit an Expression **112**
- Preview the Report **113**

1 Introduction

Access the Report Builder 5

Access Report Builder PDF Templates 6

Before You Begin 7

The Quantitative Analysis Report Builder Familiarization Guide presents step-by-step exercises to help you learn to use the Report Builder program. You can do these exercises with the demo batches DrugsofAbuse and LC-QTOF Pesticide, shipped with the system in the **Data** folder of your installation disk, or with data you acquire.

Access the Report Builder

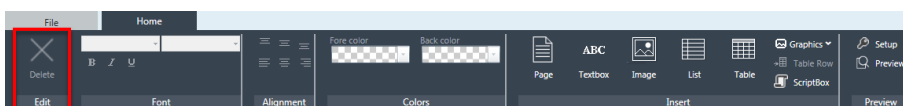
The MassHunter Quantitative Analysis Report Builder software is installed with MassHunter Quantitative Analysis.

To access the Report Builder, go to **Start > Agilent MassHunter > Report Builder**. To create a shortcut icon on your desktop, right-click **Report Builder** in the **Start** menu and select **Send to > Desktop**.

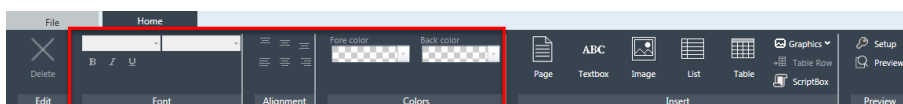
Report Builder menu

The Report Builder menu contains six sections: **Edit**, **Font**, **Alignment**, **Colors**, **Insert**, and **Preview**.

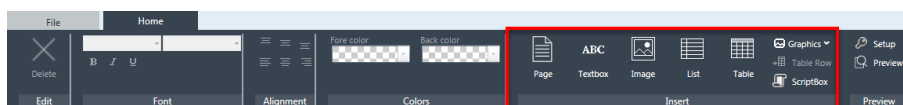
- Use the **Edit** section to **Delete** an element.



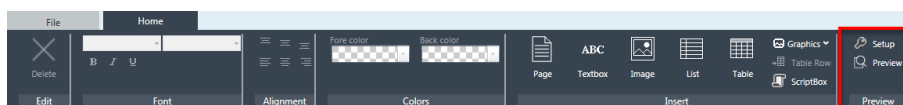
- Use the **Font**, **Alignment**, and **Colors** sections to set the font styles, results alignment, and the fore color and back color in the report template.



- Use the **Insert** section to insert **Page**, **Textbox**, **Image**, **List**, **Table**, and **Scriptbox** elements into a report template.



- Use the **Preview** section to **Setup** the Quantitative Analysis batch file that can then be used to **Preview** the report output.

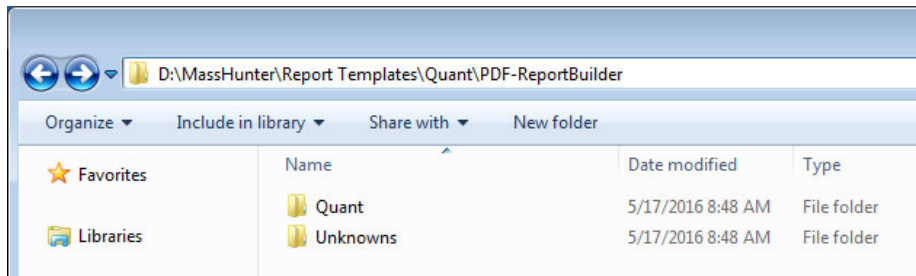


1 Introduction

Access Report Builder PDF Templates

Access Report Builder PDF Templates

Report Builder PDF Templates are located at: **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder.**



In this directory there are PDF templates for **Quant** analysis and **Unknowns** analysis. Each of these templates can be opened and edited in the Report Builder.

Example report outputs are provided on the MassHunter Quantitative Analysis Shipping DVD at **\Supplemental\MassHunter\Report Templates\Quant\ Index.**

1 Introduction

Before You Begin

Before You Begin

Either back up the PDF templates before you modify them, or immediately save the PDF template as a new template. Modified or damaged PDF templates can only be restored by re-installing the MassHunter Quantitative Analysis software.

2 Report Template Workflows

Quantitative Analysis - Calibration Report Template 9

Unknowns Analysis - Area Percent Report Template 50

In this section, you will create example report templates from scratch to give you a better understanding of how to use the MassHunter Quantitative Analysis Report Builder.

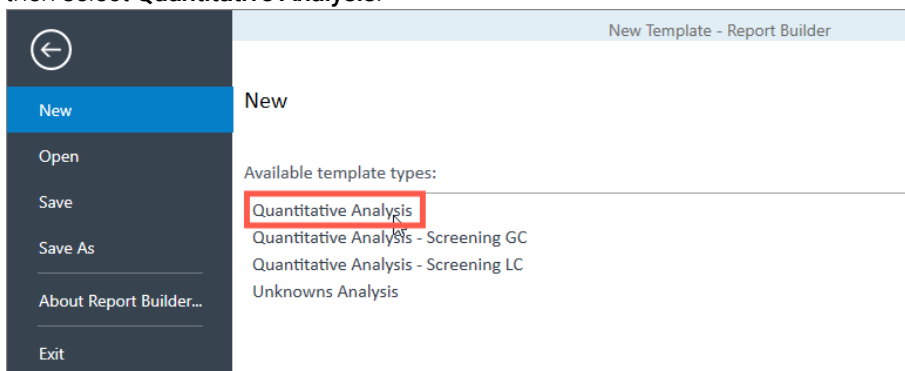
Quantitative Analysis - Calibration Report Template

For this example, you will create a Calibration Report by completing the following procedures.

- 1 **Create a new Quantitative Analysis report template** on page 9
- 2 **Edit the Batch information table** on page 14
- 3 **Add a Calibration Curve table with a graph** on page 17
- 4 **Add a Calibration Results table** on page 30
- 5 **Add a report title** on page 46
- 6 **Preview the report template** on page 48

Create a new Quantitative Analysis report template

To create a new Quantitative Analysis report template, select **File > New**, and then select **Quantitative Analysis**.



Each template type links to a different dataset. A dataset is an in-memory relational database that stores all of the information on and the relationships between samples, compounds, qualifiers, and calibration levels.

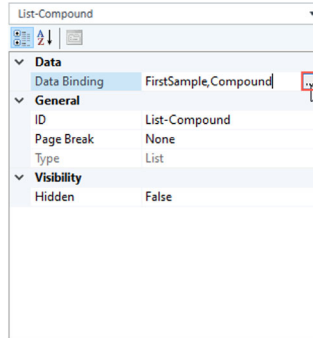
In a Quantitative Analysis dataset, a one-to-many (1:n) relationship exists between the following tables:

- Each **Batch** (sample) has one or more **TargetCompounds**.

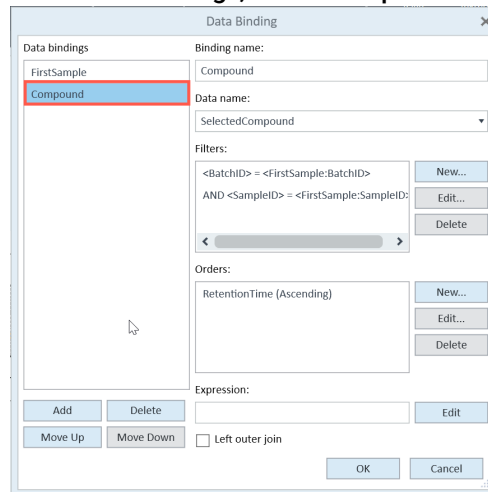
2 Report Template Workflows

Insert a list

- 3 Edit the **Data Binding** filters. Click the **Data Binding** field in the Properties pane, then click the ellipsis.



- a Under **Data bindings**, select **Compound**.



2 Report Template Workflows

Insert a list

b Under **Filters**, click **New**.

The screenshot shows the 'Data Binding' dialog box. On the left, a list of data bindings includes 'FirstSample' and 'Compound'. The 'Filters' section contains two filters: '<BatchID> = <FirstSample:BatchID>' and 'AND <SampleID> = <FirstSample:SampleID>'. A red box highlights the 'New...' button next to the first filter. Below the filters are 'Orders' and 'Expression' sections, each with 'New...', 'Edit...', and 'Delete' buttons. At the bottom, there are 'Add', 'Delete', 'Move Up', 'Move Down', and 'Left outer join' options, along with 'OK' and 'Cancel' buttons.

NOTE

An ID is a unique identifier for a table row in a Dataset. IDs such as BatchID, SampleID, CompoundID, etc. have system-assigned numbers. The combination of IDs can rapidly locate and link to the unique table row in a Quantitative Analysis dataset.

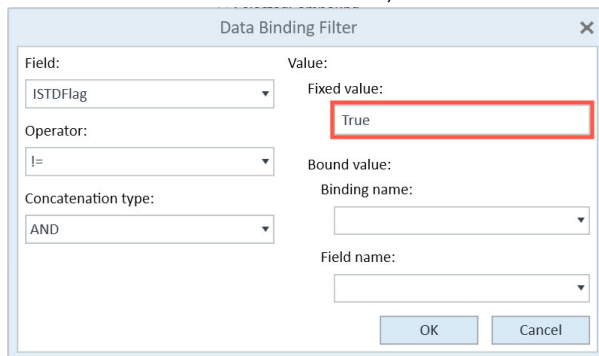
c Select **ISTDFlag** from the **Field** list, and select **!=** from the **Operator** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'ISTDFlag' and the 'Operator' dropdown is set to '!=', both highlighted with red boxes. The 'Concatenation type' is set to 'AND'. The 'Value' section has empty fields for 'Fixed value', 'Bound value', and 'Field name'. 'OK' and 'Cancel' buttons are at the bottom.

2 Report Template Workflows

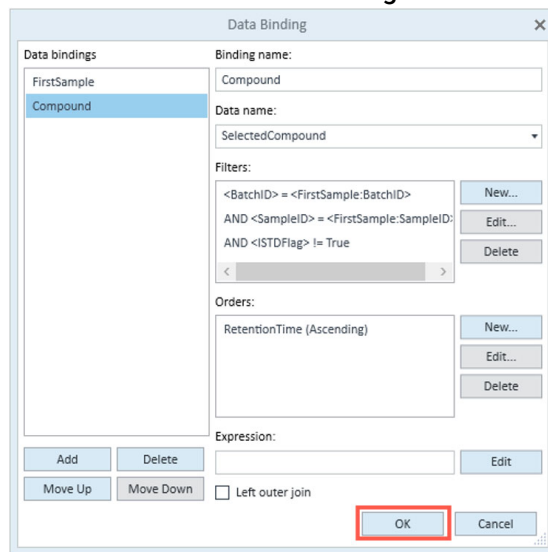
Insert a list

d Enter **True** in the **Fixed value** field, and click **OK**.



The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'ISTDFlag'. The 'Fixed value' field contains the text 'True', which is highlighted with a red rectangular box. The 'Operator' dropdown is set to '!='. The 'Concatenation type' dropdown is set to 'AND'. The 'Value', 'Bound value', 'Binding name', and 'Field name' fields are empty. The 'OK' and 'Cancel' buttons are at the bottom right.

e Click **OK** to close the **Data Binding** window.

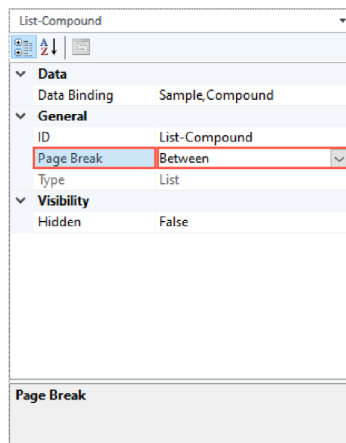


The screenshot shows the 'Data Binding' dialog box. The 'Data bindings' list on the left contains 'FirstSample' and 'Compound', with 'Compound' selected. The 'Binding name' is 'Compound'. The 'Data name' dropdown is set to 'SelectedCompound'. The 'Filters' section contains three filter expressions: '<BatchID> = <FirstSample:BatchID>', 'AND <SampleID> = <FirstSample:SampleID>', and 'AND <ISTDFlag> != True'. The 'Orders' section contains 'RetentionTime (Ascending)'. The 'Expression' field is empty. The 'Left outer join' checkbox is unchecked. The 'OK' button is highlighted with a red rectangular box. The 'Add', 'Delete', 'Move Up', and 'Move Down' buttons are on the left, and 'New...', 'Edit...', and 'Delete' buttons are on the right.

2 Report Template Workflows

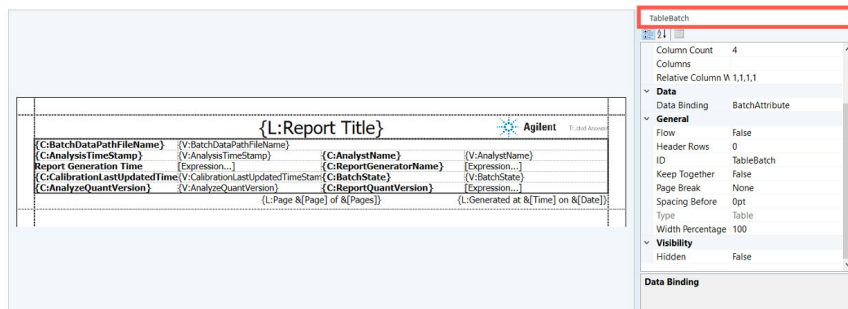
Edit the Batch information table

- 4 In the Properties pane, under General, click **Page Break**, then select **Between** from the list.



Edit the Batch information table

- 1 Select **TableBatch** from the drop-down menu at the top of the Properties pane. This highlights the table filled with batch information in your report. Each cell contains a caption, value, or expression.



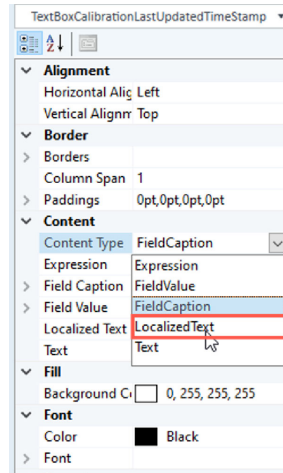
A table contains a header row, rows, and cells. Table cells can include the following:

- **(C)** - Caption (same term that is displayed in the Quantitative Analysis software)
- **(L)** - Localized text
- **(V)** - Value (same number that is displayed in the Quantitative Analysis software)
- **(E)** - Expressions (can be used to format numbers and data as desired)

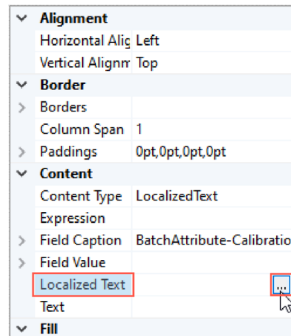
2 Report Template Workflows

Edit the Batch information table

- The caption **{C:CalibrationLastUpdatedTimeStamp}** is too long to fit in its cell. Select it, click **Content Type** in the Properties pane, and select **LocalizedText** from the drop-down menu.



- Click **Localized Text**, then click the ellipsis.

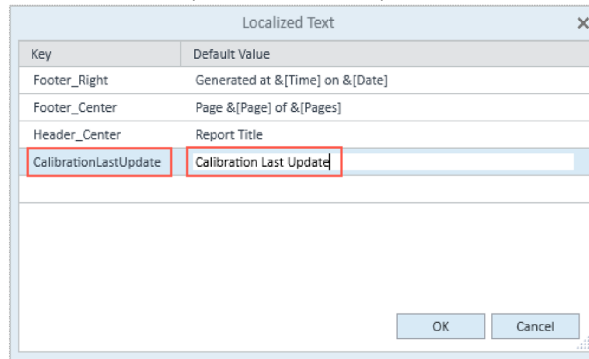


- On the bottom row, enter **CalibrationLastUpdate** in the **Key** column, and **Calibration Last Update** in the **Default Value** column.

2 Report Template Workflows

Edit the Batch information table

- 5 Click **OK**. The caption has been updated in the table.

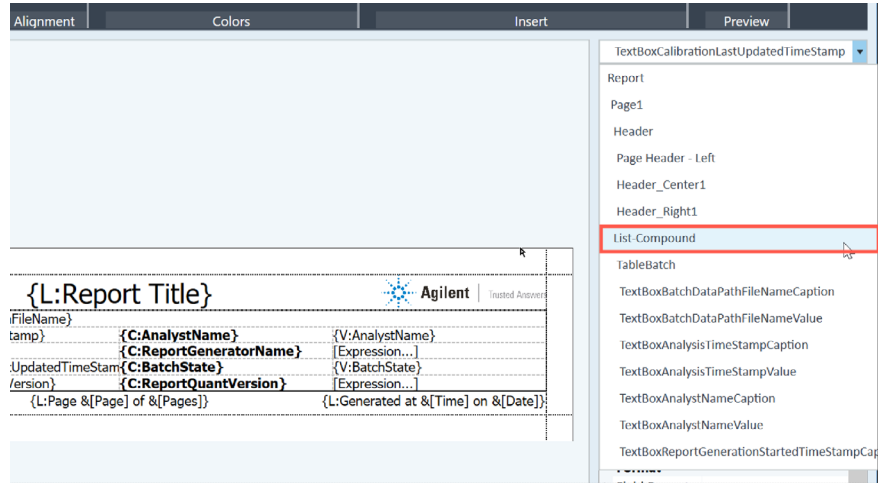


2 Report Template Workflows

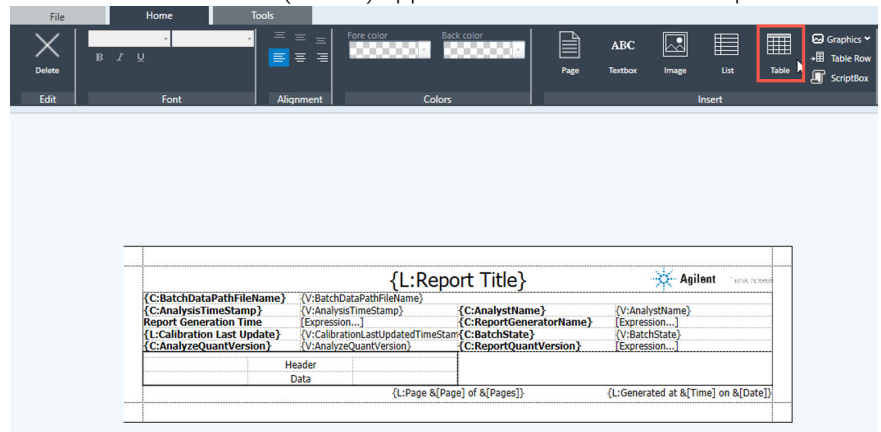
Add a Calibration Curve table with a graph

Add a Calibration Curve table with a graph

- 1 Select **List-Compound** from the drop-down menu at the top of the Properties pane.



- 2 Click **Table**. A new table (Table1) appears at the bottom of the report.



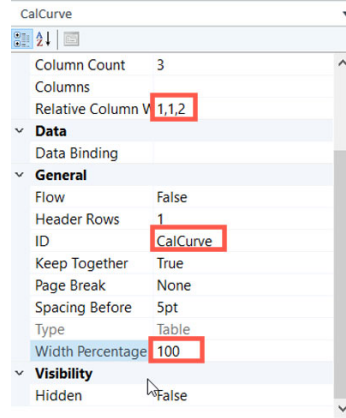
In the properties pane, make the following changes:

- Change the **Relative Column Width** to 1,1,2 to resize the columns in relation to the table width
- Enter CalCurve in the **ID** field to update the table name

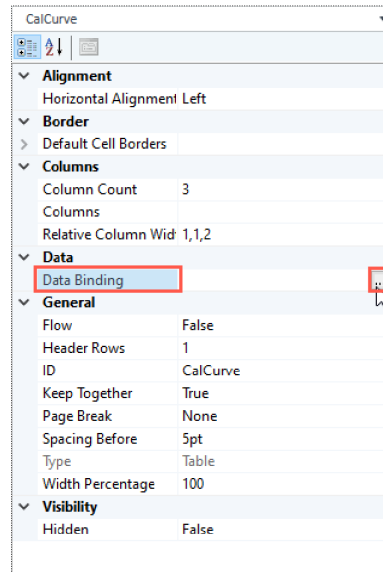
2 Report Template Workflows

Add a Calibration Curve table with a graph

- Enter 100 in the **Width Percentage** field to resize the table width



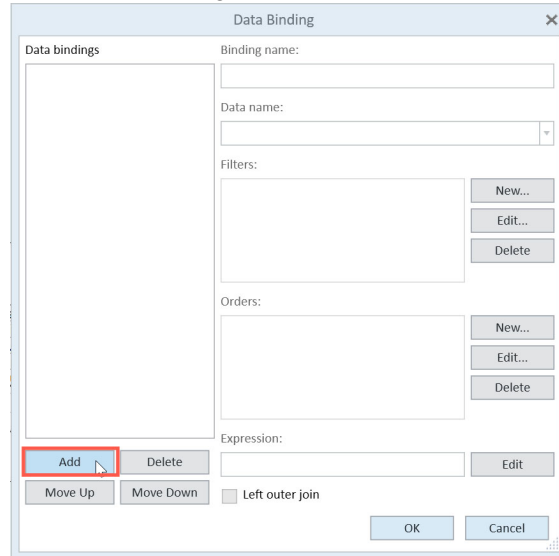
- 3 To add binding data to the table, click **Data Binding** in the Properties pane, then click the ellipsis.



2 Report Template Workflows

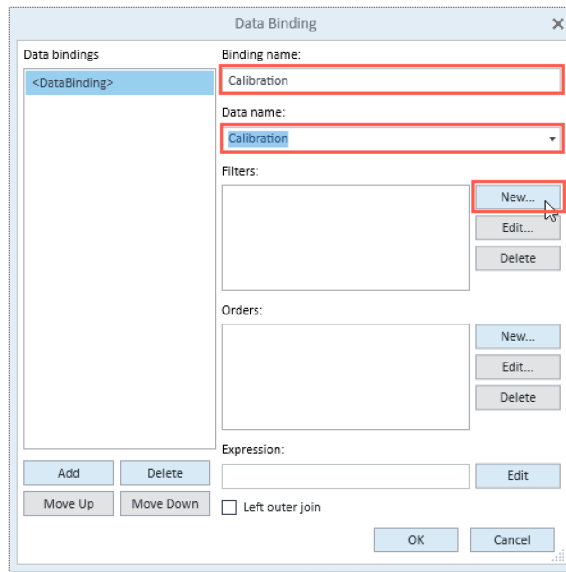
Add a Calibration Curve table with a graph

4 Under **Data Bindings**, click **Add**.



The screenshot shows the 'Data Binding' dialog box. On the left, there is a list titled 'Data bindings' which is currently empty. Below this list are four buttons: 'Add', 'Delete', 'Move Up', and 'Move Down'. The 'Add' button is highlighted with a red rectangle. To the right of the list, there are several input fields: 'Binding name:' (empty), 'Data name:' (empty), 'Filters:' (empty), 'Orders:' (empty), and 'Expression:' (empty). Each of these fields has a corresponding 'New...', 'Edit...', or 'Delete' button. At the bottom right, there are 'OK' and 'Cancel' buttons.

5 Enter **Calibration** for the **Binding name** and select **Calibration** from the **Data name** list. In the **Filters** section, click **New**.



The screenshot shows the 'Data Binding' dialog box after the first step. The 'Data bindings' list now contains one entry: '<DataBinding>'. The 'Binding name:' field contains the text 'Calibration'. The 'Data name:' dropdown menu is open, showing 'Calibration' selected. In the 'Filters:' section, the 'New...' button is highlighted with a red rectangle. The 'Orders:' section is empty. The 'Expression:' field is empty. The 'Left outer join' checkbox is unchecked. The 'Add', 'Delete', 'Move Up', and 'Move Down' buttons are now disabled. The 'New...', 'Edit...', and 'Delete' buttons for the 'Filters' and 'Orders' sections are also disabled. The 'OK' and 'Cancel' buttons are at the bottom right.

2 Report Template Workflows

Add a Calibration Curve table with a graph

- 6 Select **SampleID** from the **Field** list and **=** from the **Operator** list. Select **Compound** from the **Binding name** list, and select **SampleID** from the **Field name** list. Click **OK**.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'SampleID', the 'Operator' dropdown is set to '=', the 'Binding name' dropdown is set to 'Compound', and the 'Field name' dropdown is set to 'SampleID'. The 'Fixed value' and 'Concatenation type' fields are empty. The 'OK' button is highlighted with a mouse cursor.

- 7 Add a data binding filter for each of the following configurations:
 - a Select **CompoundID** from the **Field** list, select **=** from the **Operator** list, select **AND** from the **Concatenation type** list, select **Compound** from the **Binding name** list, and select **CompoundID** from the **Field name** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'CompoundID', the 'Operator' dropdown is set to '=', the 'Concatenation type' dropdown is set to 'AND', the 'Binding name' dropdown is set to 'Compound', and the 'Field name' dropdown is set to 'CompoundID'. The 'Fixed value' field is empty. The 'OK' and 'Cancel' buttons are visible at the bottom.

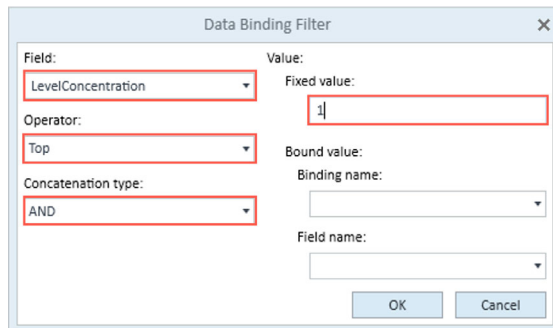
- b Select **CalibrationType** from the **Field** list, select **!=** from the **Operator** list, select **AND** from the **Concatenation type** list, and select **Method** from the **Fixed value** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'CalibrationType', the 'Operator' dropdown is set to '!=', the 'Concatenation type' dropdown is set to 'AND', and the 'Fixed value' text box contains 'Method'. The 'Binding name' and 'Field name' dropdowns are empty. The 'OK' and 'Cancel' buttons are visible at the bottom.

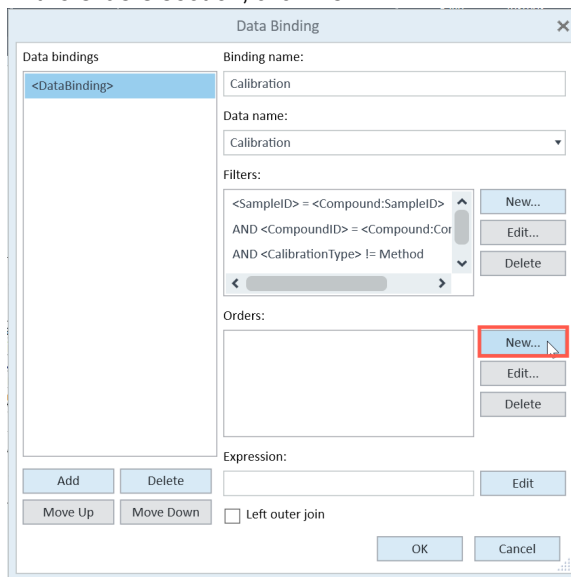
2 Report Template Workflows

Add a Calibration Curve table with a graph

- c Select **LevelConcentration** from the **Field** list, select **Top** from the **Operator** list, select **AND** from the **Concatenation type** list, and enter **1** for the **Fixed value**.



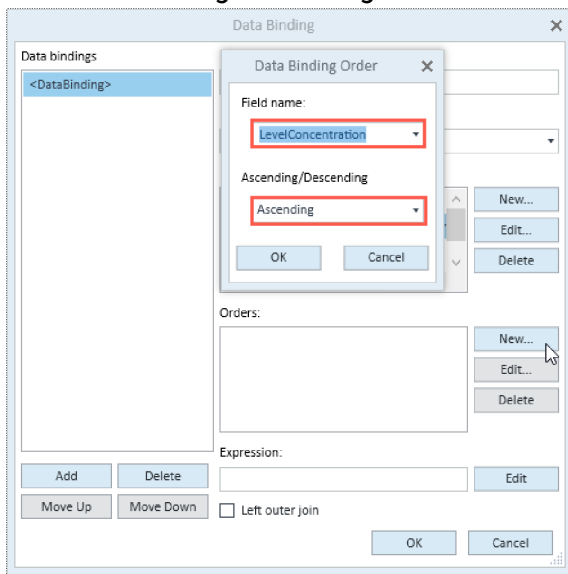
- 8 In the **Orders** section, click **New**.



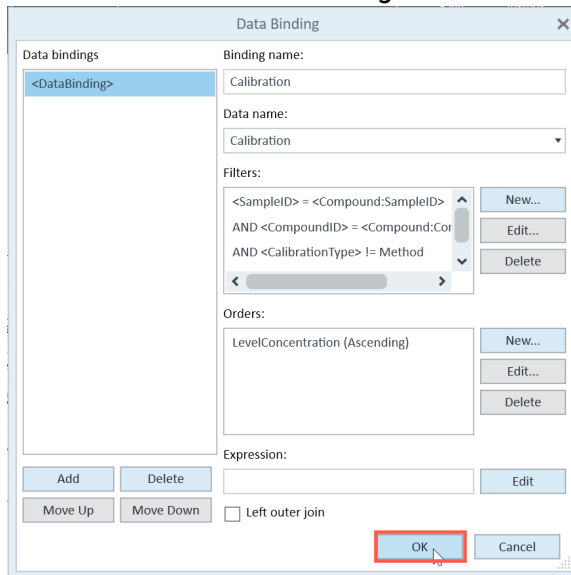
2 Report Template Workflows

Add a Calibration Curve table with a graph

- 9 Select **LevelConcentration** from the **Field name** list, and select **Ascending** from the **Ascending/Descending** list. Click **OK**.



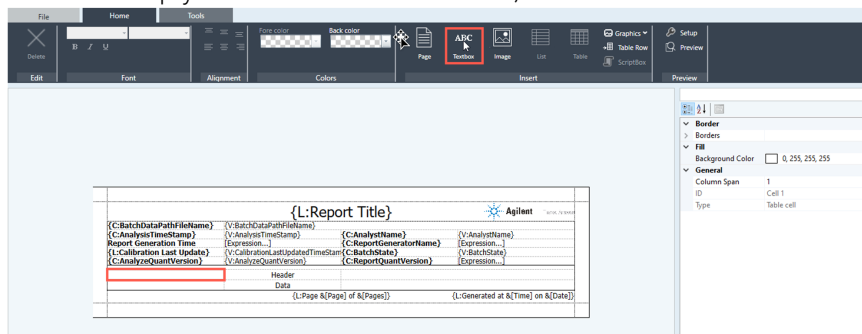
- 10 Click **OK** to close the **Data Binding** window.



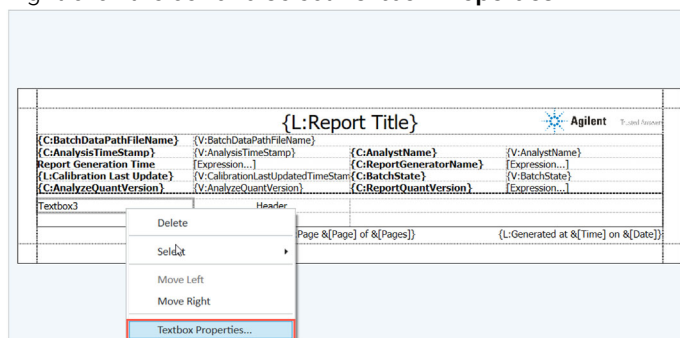
2 Report Template Workflows

Add a Calibration Curve table with a graph

11 Select the empty cell to the left of the **Header**, and click **Textbox**.



12 Right-click the cell and select **Textbox Properties**.



2 Report Template Workflows

Add a Calibration Curve table with a graph

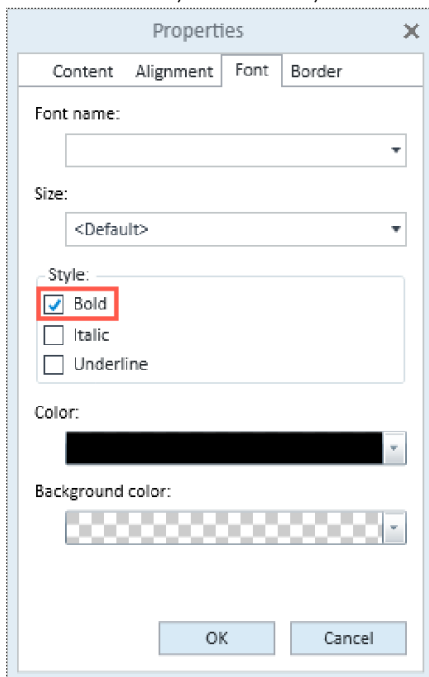
- 13 Select **Field value**, select **Compound** from the **Binding name** list, and then select **CompoundName** from the **Field** list.

The screenshot shows a 'Properties' dialog box with a close button (X) in the top right corner. The dialog has four tabs: 'Content', 'Alignment', 'Font', and 'Border'. The 'Content' tab is selected. There are four radio button options: 'Localized text', 'Text', 'Field value', and 'Field caption'. The 'Field value' option is selected and highlighted with a red box. Below the radio buttons, there are two dropdown menus. The first dropdown is labeled 'Binding name:' and has 'Compound' selected, also highlighted with a red box. The second dropdown is labeled 'Field:' and has 'CompoundName' selected, also highlighted with a red box. Below these are two more dropdown menus for 'Field caption', both of which are empty. At the bottom of the dialog are 'OK' and 'Cancel' buttons.

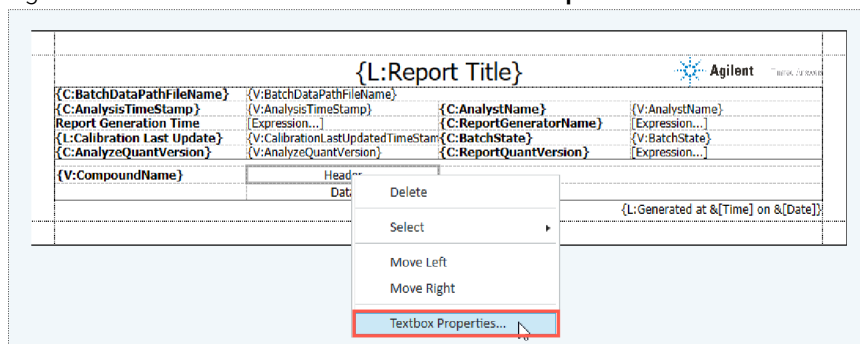
2 Report Template Workflows

Add a Calibration Curve table with a graph

14 On the **Font** tab, select **Bold**, and click **OK**.



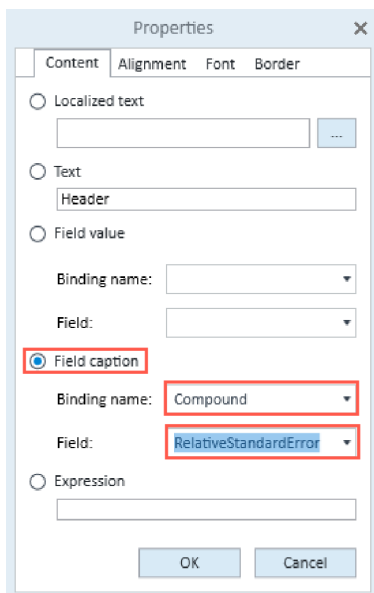
15 Right-click the **Header** cell and select **Textbox Properties**.



2 Report Template Workflows

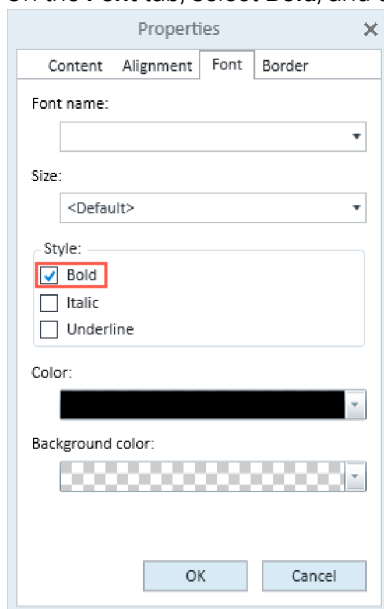
Add a Calibration Curve table with a graph

- 16 Select **Field Caption**, select **Compound** from the **Binding name** list, and select **RelativeStandardError** from the **Field** list.



The screenshot shows the 'Properties' dialog box with the 'Field caption' tab selected. The 'Field caption' radio button is highlighted with a red box. The 'Binding name' dropdown is set to 'Compound' and the 'Field' dropdown is set to 'RelativeStandardError', both also highlighted with red boxes. The 'OK' and 'Cancel' buttons are at the bottom.

- 17 On the **Font** tab, select **Bold**, and click **OK**.

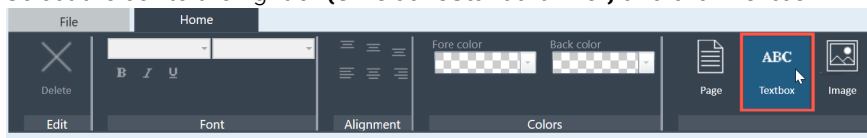


The screenshot shows the 'Properties' dialog box with the 'Font' tab selected. The 'Bold' checkbox in the 'Style' section is checked and highlighted with a red box. The 'OK' and 'Cancel' buttons are at the bottom.

2 Report Template Workflows

Add a Calibration Curve table with a graph

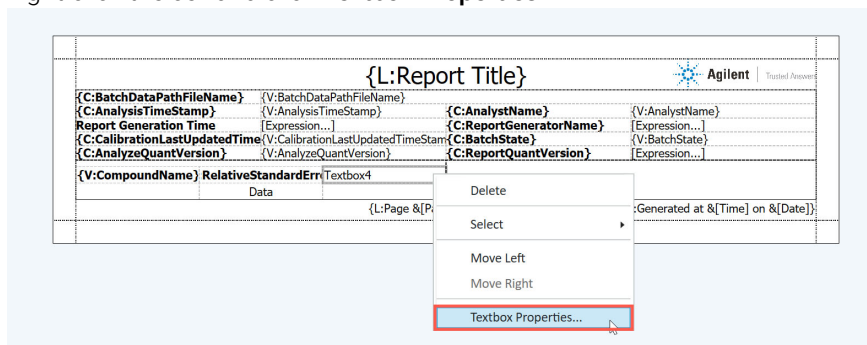
18 Select the cell to the right of **{C:RelativeStandardError}** and click **Textbox**.



The screenshot shows the Microsoft Word ribbon with the 'Home' tab selected. The 'Page' group contains the 'Page', 'Textbox', and 'Image' buttons. The 'Textbox' button is highlighted with a red box. Below the ribbon, a portion of a report template is visible, showing a table with various fields and a red box around a cell.

		{L:Report Title}
{C:BatchDataPathFileName}	{V:BatchDataPathFileName}	
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:AnalystName}
Report Generation Time	[Expression...]	{C:ReportGeneratorName}
{C:CalibrationLastUpdatedTime}	{V:CalibrationLastUpdatedTime}	{C:BatchState}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}	{C:ReportQuantVersion}
{V:CompoundName}	RelativeStandardErr	
	Data	
		{L:Page &[Page] of &[Pages]}

19 Right-click the cell and click **Textbox Properties**.



The screenshot shows the same report template as in the previous image. A right-click context menu is open over the cell containing the text 'RelativeStandardErr'. The 'Textbox Properties...' option is highlighted with a red box.

		{L:Report Title}	Agilent	Trusted Access
{C:BatchDataPathFileName}	{V:BatchDataPathFileName}		{C:AnalystName}	{V:AnalystName}
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}		{C:ReportGeneratorName}	[Expression...]
Report Generation Time	[Expression...]		{C:BatchState}	{V:BatchState}
{C:CalibrationLastUpdatedTime}	{V:CalibrationLastUpdatedTime}		{C:ReportQuantVersion}	[Expression...]
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}			
{V:CompoundName}	RelativeStandardErr	Textbox4		
	Data			
		{L:Page &[Page] of &[Pages]}		Generated at &[Time] on &[Date]

2 Report Template Workflows

Add a Calibration Curve table with a graph

- 20 Select **Field Value**, select **Compound** from the **Binding name** list, and **RelativeStandardError** from the **Field** list. Click **OK**.

The screenshot shows the 'Properties' dialog box for 'Textbox4'. The 'Field value' radio button is selected and highlighted with a red box. The 'Binding name' dropdown menu is set to 'Compound' and the 'Field' dropdown menu is set to 'RelativeStandardError', both also highlighted with red boxes. Other options like 'Localized text', 'Text', 'Field caption', and 'Expression' are unselected. The 'OK' and 'Cancel' buttons are at the bottom.

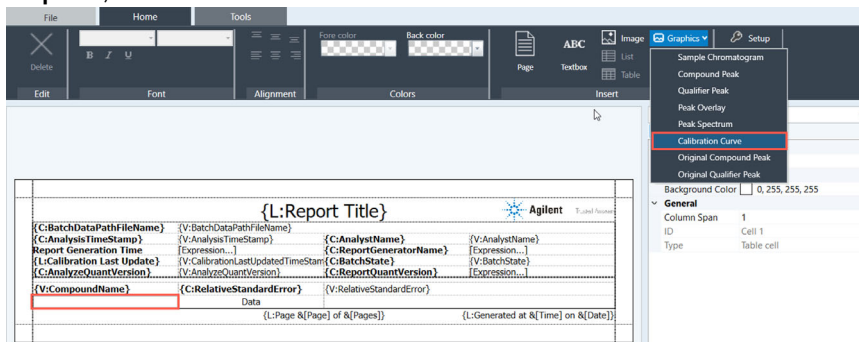
- 21 To use the same format in the current cell as is found in the Quantitative Analysis software, double-click **Field Format** in the **Properties** pane, and select **Compound** from the **Binding Name** list, and **RelativeStandardError** from the **Field Name** list.

The screenshot shows the 'Properties' pane for 'Textbox4'. The 'Field Format' section is expanded and highlighted with a red box. Within this section, the 'Binding Name' dropdown is set to 'Compound' and the 'Field Name' dropdown is set to 'RelativeStandardError', both highlighted with red boxes. Other sections like 'Content', 'Fill', 'Font', and 'General' are also visible.

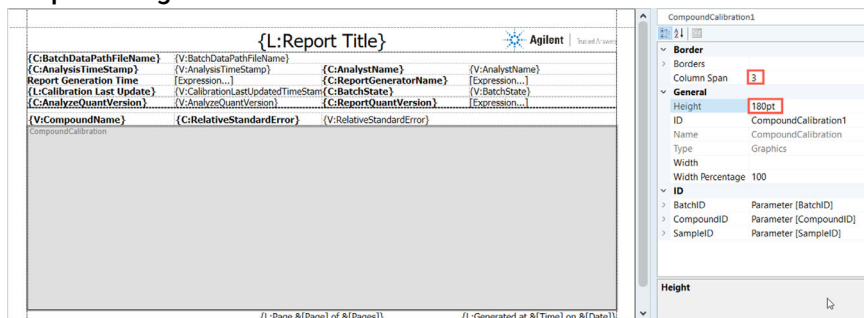
2 Report Template Workflows

Add a Calibration Curve table with a graph

- 22 To add a graphic, select the empty cell below **{V:CompoundName}**, click **Graphics**, and select **Calibration Curve**.



- 23 To resize the graphic, in the Properties pane, enter **3** for **Column Span** and **180pt** for **Height**.



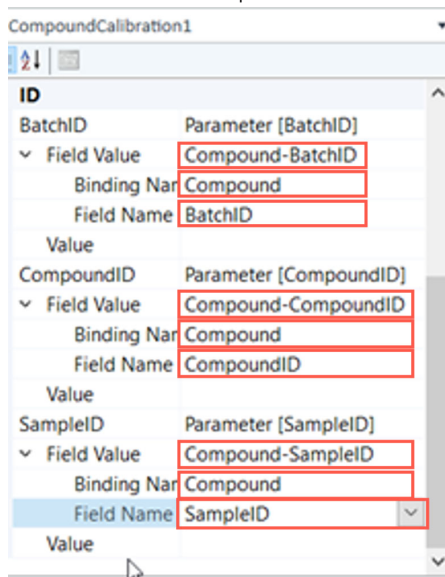
ID fields like BatchID, SampleID, and CompoundID have system-assigned numbers which help to rapidly locate and link to unique table rows in a Quantitative Analysis dataset. To edit the ID fields, expand the ID section in the Properties pane. Enter the values for each ID:

- BatchID
 - **Field Value:** Compound-BatchID
 - **Binding Name:** Compound
 - **Field Name:** BatchID
- CompoundID
 - **Field Value:** Compound-CompoundID
 - **Binding Name:** Compound
 - **Field Name:** CompoundID
- SampleID
 - **Field Value:** Compound-SampleID

2 Report Template Workflows

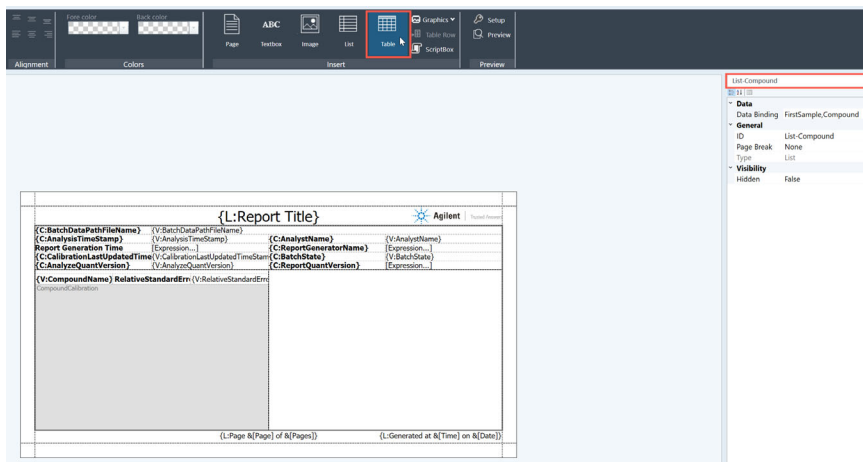
Add a Calibration Results table

- **Binding Name:** Compound
- **Field Name:** SampleID



Add a Calibration Results table

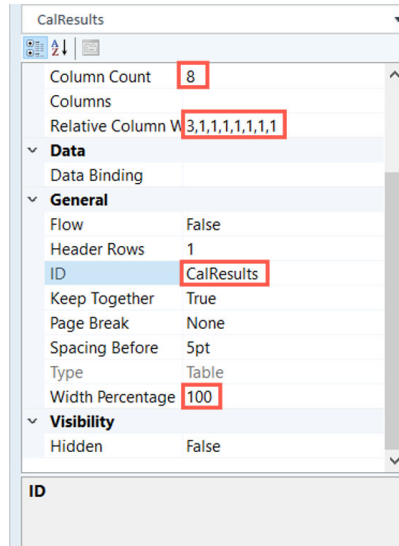
- 1 Select **List-Compound** from the list at the top of the Properties pane, and click **Table**.



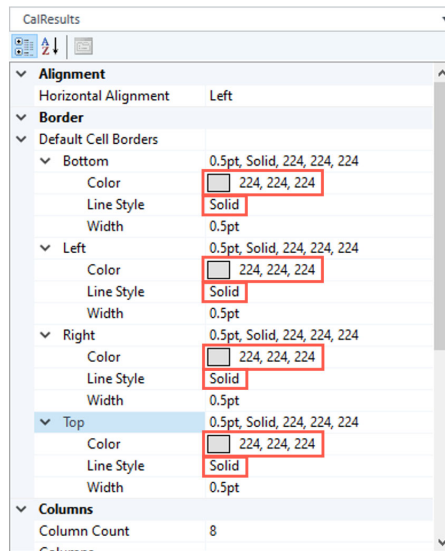
2 Report Template Workflows

Add a Calibration Results table

- In the Properties pane, adjust the number of columns by changing the **Column Count** to **8**, and adjust the widths of the columns by changing **Relative Column Widths** to **3,1,1,1,1,1,1,1**. Enter **CalResults** for **ID**, and **100** for **Width Percentage**.



- To add a border to the table, select **Solid** from the **Line Style** list and select a **Color** for each of the cell borders.

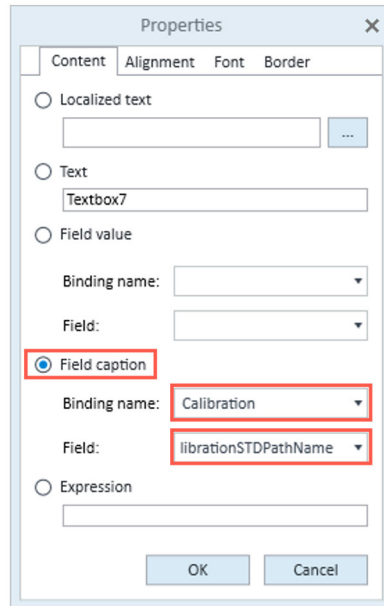


- Repeat steps 3-10 from the section **Add a Calibration Curve table with a graph** on page 17 to add binding data to the table.

2 Report Template Workflows

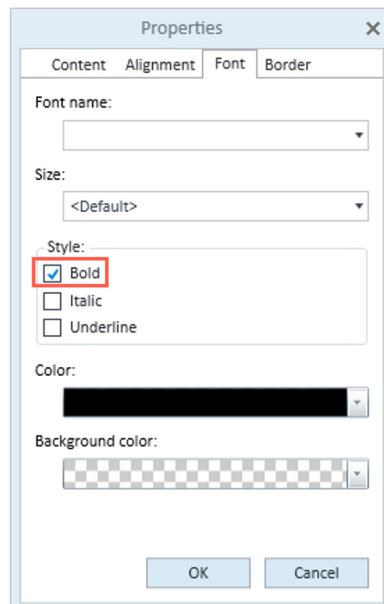
Add a Calibration Results table

- 6 Right-click the cell and select **Textbox Properties**. Select **Field Caption**, select **Calibration** from the **Binding name** list, and **CalibrationSTDPATHName** from the **Field** list.



The screenshot shows the 'Properties' dialog box with the 'Field Caption' tab selected. The 'Field caption' radio button is selected and highlighted with a red box. The 'Binding name' dropdown is set to 'Calibration' and the 'Field' dropdown is set to 'librationSTDPATHName', both also highlighted with red boxes. The 'Content' tab is active, and the 'Text' option is selected. The 'Field value' section is empty. The 'OK' and 'Cancel' buttons are at the bottom.

- 7 On the **Font** tab, select **Bold**. Click **OK**.



The screenshot shows the 'Properties' dialog box with the 'Font' tab selected. The 'Style' section has the 'Bold' checkbox checked and highlighted with a red box. The 'Font name' dropdown is empty, 'Size' is '<Default>', 'Color' is black, and 'Background color' is a checkerboard pattern. The 'Content' tab is active. The 'OK' and 'Cancel' buttons are at the bottom.

2 Report Template Workflows

Add a Calibration Results table

- 8 Select the first cell in the second row of the table, and click **Textbox**.

The screenshot shows the Report Builder interface. At the top, there is a ribbon with various options. The 'Insert' tab is active, and the 'Textbox' option is highlighted with a red box. Below the ribbon, a report template is displayed. The template includes a header section with the following fields:

{L:Report Title}		Agilent Trusted Answers	
{C:BatchDataPathFileName}	{V:BatchDataPathFileName}	{C:AnalystName}	{V:AnalystName}
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:ReportGeneratorName}	{V:ReportGeneratorName}
Report Generation Time	[Expression...]	{C:BatchState}	{V:BatchState}
{C:CalibrationLastUpdatedTime}	{V:CalibrationLastUpdatedTime}	{C:ReportQuantVersion}	{V:ReportQuantVersion}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}		

Below the header is a table with the following structure:

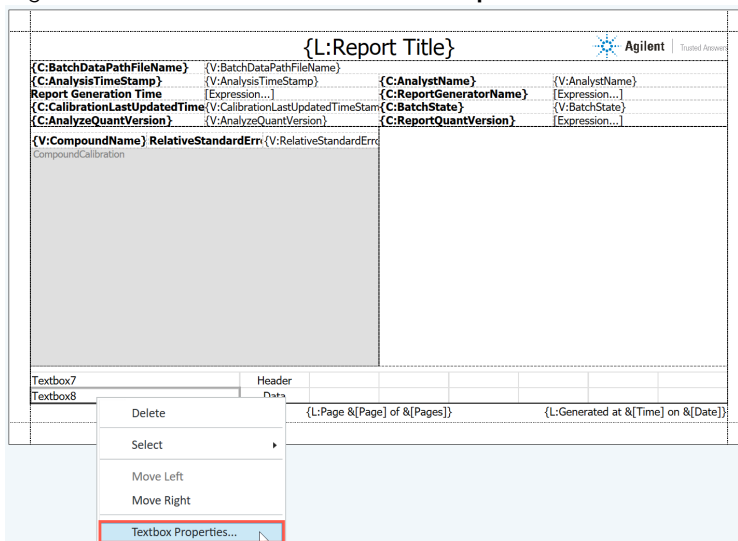
{V:CompoundName}	RelativeStandardErrr	{V:RelativeStandardErrr}
CompoundCalibration		

At the bottom of the table, the first cell of the second row is highlighted with a red box and labeled 'Textbox7'. The table has a 'Header' row and a 'Data' row. The footer of the report includes the page number and generation time: {L:Page &[Page] of &[Pages]} and {L:Generated at &[Time] on &[Date]}.

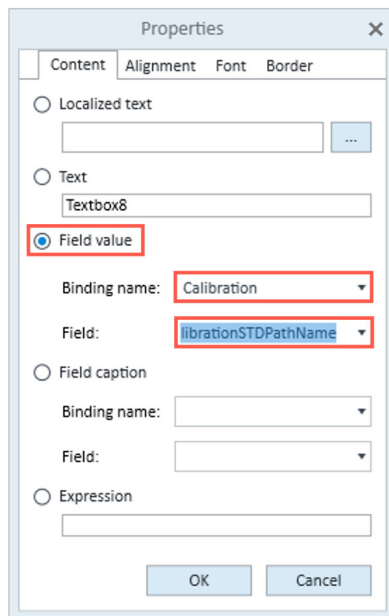
2 Report Template Workflows

Add a Calibration Results table

9 Right-click the cell and select **Textbox Properties**.



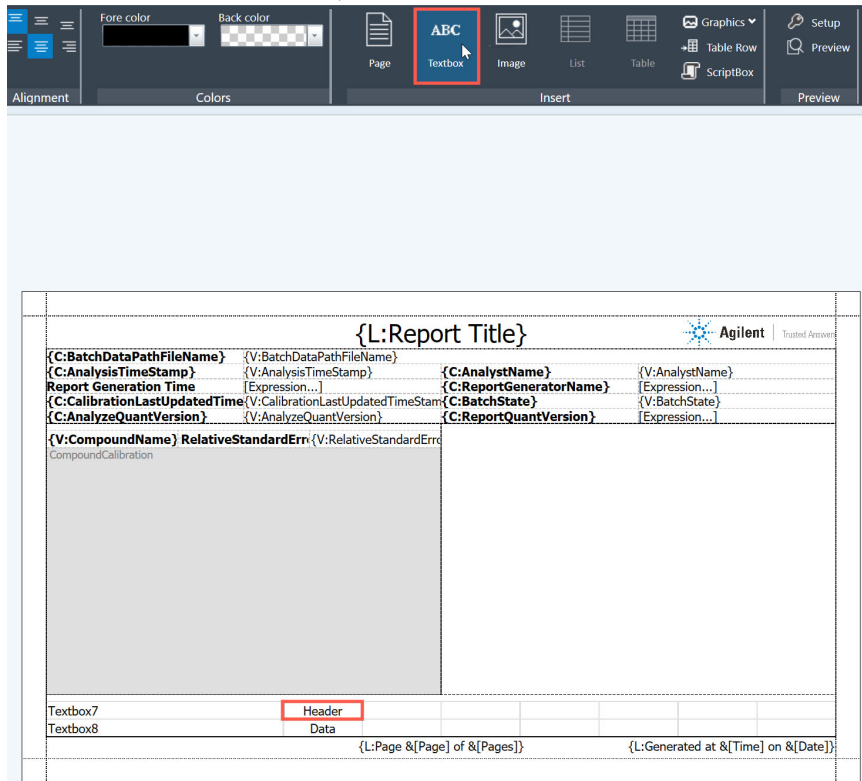
10 Select **Field Value**, select **Calibration** from the **Binding name** list, and select **CalibrationSTDPATHName** from the **Field** list. Click **OK**.



2 Report Template Workflows

Add a Calibration Results table

11 Select the cell labeled **Header**, and click **Textbox**.

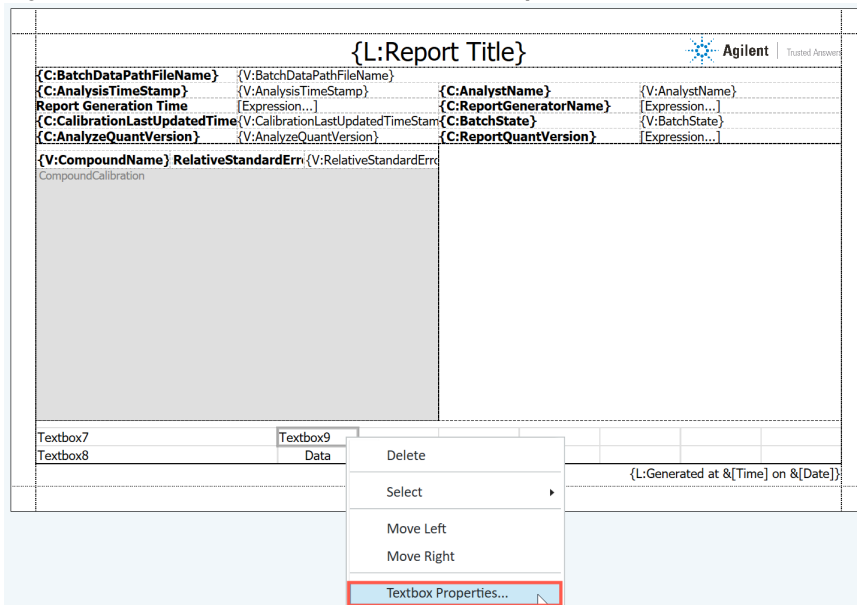


The screenshot displays the Report Builder software interface. At the top, a ribbon menu is visible with sections for 'Alignment', 'Colors', 'Insert', and 'Preview'. The 'Insert' section contains icons for 'Page', 'Textbox', 'Image', 'List', and 'Table'. The 'Textbox' icon, which shows the letters 'ABC', is highlighted with a red rectangular box. Below the ribbon, a report template is shown. The template includes a header area with the text '{L:Report Title}' and the Agilent logo. Below the header is a table with several rows of metadata, including fields like '{C:BatchDataPathFileName}', '{C:AnalysisTimeStamp}', and '{C:ReportGeneratorName}'. A large table with a grey background is also present, with the header row containing '{V:CompoundName}' and 'RelativeStandardError'. At the bottom of the report, there are two textboxes labeled 'Textbox7' and 'Textbox8'. 'Textbox7' contains the word 'Header' and is highlighted with a red box. 'Textbox8' contains the word 'Data'. The footer of the report includes page and date information: '{L:Page &[Page] of &[Pages]}' and '{L:Generated at &[Time] on &[Date]}'. The 'Agilent' logo and 'Trusted Advisor' text are visible in the top right corner of the report area.

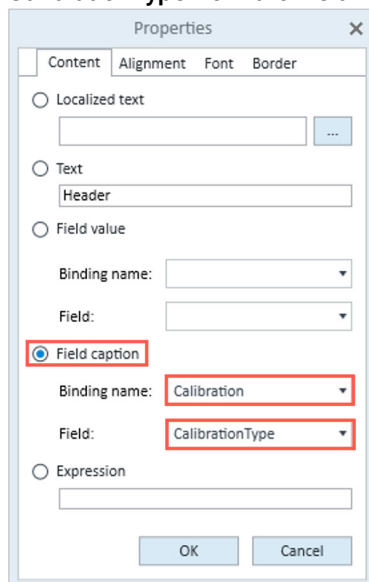
2 Report Template Workflows

Add a Calibration Results table

12 Right-click the textbox, and select **Textbox Properties**.



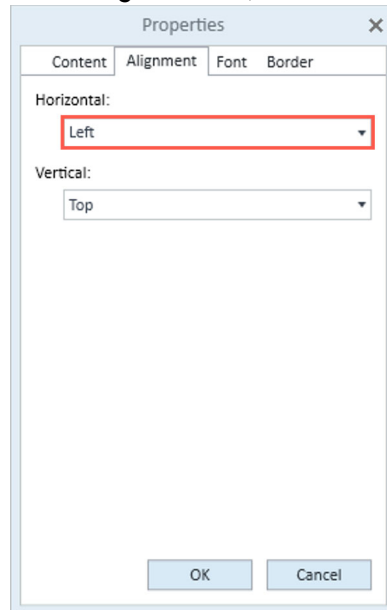
13 Select **Field Caption**, select **Calibration** from the **Binding name** list, and **CalibrationType** from the **Field** list.



2 Report Template Workflows

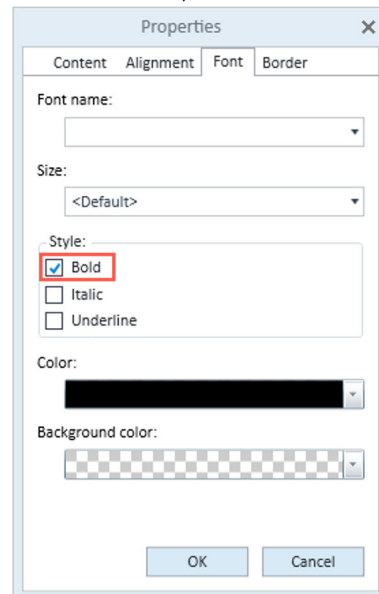
Add a Calibration Results table

14 On the **Alignment** tab, select **Left** from the **Horizontal** list.



The screenshot shows the 'Properties' dialog box with the 'Alignment' tab selected. The 'Horizontal' dropdown menu is set to 'Left' and is highlighted with a red rectangular box. The 'Vertical' dropdown menu is set to 'Top'. At the bottom of the dialog are 'OK' and 'Cancel' buttons.

15 On the **Font** tab, select **Bold**. Click **OK**.



The screenshot shows the 'Properties' dialog box with the 'Font' tab selected. The 'Style' section has the 'Bold' checkbox checked and highlighted with a red rectangular box. The 'Font name' dropdown is empty, 'Size' is '<Default>', 'Color' is black, and 'Background color' is a checkerboard pattern. At the bottom of the dialog are 'OK' and 'Cancel' buttons.

2 Report Template Workflows

Add a Calibration Results table

16 Select the cell labeled **Data**, and click **Textbox**.

The screenshot shows the Report Builder interface. At the top, the 'Insert' menu is open, and the 'Textbox' option is highlighted with a red box. Below the menu, a report template is displayed. The template includes a header section with the title '{L:Report Title}' and the Agilent logo. Below the header, there is a table with the following content:

{C:BatchDataPathFileName}	{V:BatchDataPathFileName}	{C:AnalystName}	{V:AnalystName}
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:ReportGeneratorName}	{V:ReportGeneratorName}
Report Generation Time	[Expression...]	{C:BatchState}	{V:BatchState}
{C:CalibrationLastUpdatedTime}	{V:CalibrationLastUpdatedTime}	{C:ReportQuantVersion}	{V:ReportQuantVersion}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}		

Below the table, there is a large greyed-out area labeled 'CompoundCalibration'. At the bottom of the report, there is a table with the following content:

{C:CalibrationSTDPathName}	{C:Calibratic		
{V:CalibrationSTDPathName}	Data		

The 'Data' cell in the bottom table is highlighted with a red box. At the bottom right of the report, there is a footer with the text '{L:Page &[Page] of &[Pages]}' and '{L:Generated at &[Time] on &[Date]}'. The Agilent logo and 'Trusted Account' are visible in the top right corner of the report area.

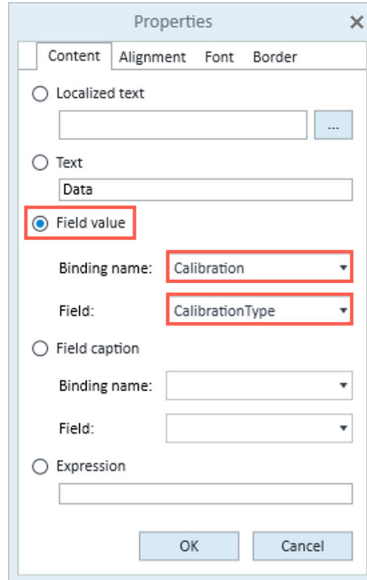
17 Right-click the cell, and select **Textbox Properties**.

The screenshot shows the Report Builder interface with the 'Textbox Properties' context menu open over the 'Data' cell. The menu options are: Delete, Select, Move Left, Move Right, and Textbox Properties... (highlighted with a red box). The report template content is the same as in the previous screenshot.

2 Report Template Workflows

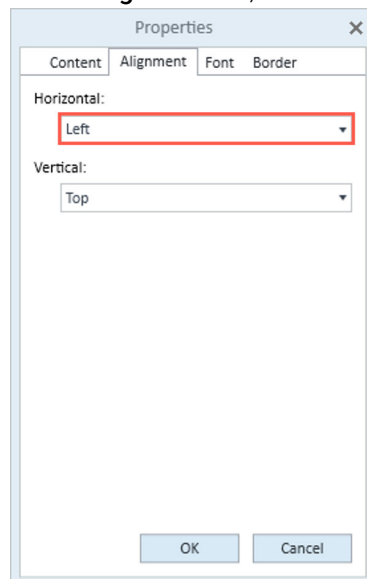
Add a Calibration Results table

- 18 Select **Field Value**, select **Calibration** from the **Binding name** list, and select **CalibrationType** from the **Field** list.



The screenshot shows the 'Properties' dialog box with the 'Field value' radio button selected. The 'Binding name' dropdown is set to 'Calibration' and the 'Field' dropdown is set to 'CalibrationType'. Both dropdowns are highlighted with a red box. The 'Field caption' and 'Expression' options are not selected. The 'Content' tab is active, and the 'Text' option is selected. The 'Data' field is empty. The 'OK' and 'Cancel' buttons are at the bottom.

- 19 On the **Alignment** tab, select **Left** from the **Horizontal** list. Click **OK**.



The screenshot shows the 'Properties' dialog box with the 'Alignment' tab selected. The 'Horizontal' dropdown is set to 'Left' and the 'Vertical' dropdown is set to 'Top'. Both dropdowns are highlighted with a red box. The 'OK' and 'Cancel' buttons are at the bottom.

2 Report Template Workflows

Add a Calibration Results table

- 20 For the cells in the third column, use the instructions in the previous steps to select **Calibration** and **LevelName** for the **Field Caption** of the cell in the first row, and **Calibration** and **LevelName** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Center**. Click **OK**.

{L:Report Title}			
{C:BatchDataPathFileName}	{V:BatchDataPathFileName}		
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:AnalystName}	{V:AnalystName}
Report Generation Time	{Expression...}	{C:ReportGeneratorName}	{Expression...}
{L:Calibration Last Update}	{V:CalibrationLastUpdatedTimeStamp}	{C:BatchState}	{V:BatchState}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}	{C:ReportQuantVersion}	{Expression...}
{V:CompoundName}	{C:RelativeStandardError}	{V:RelativeStandardError}	
CompoundCalibration			
{C:CalibrationSTDPPathName}	{C:CalibratiC:LevelName}		
{V:CalibrationSTDPPathName}	{V:CalibrationV:LevelName}		
		{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

- 21 For the cells in the fourth column, select **Calibration** and **LevelEnable** for the **Field Caption** of the cell in the first row, and **Calibration** and **LevelEnable** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and set the horizontal alignment of both cells to **Center**. Click **OK**.

{L:Report Title}			
{C:BatchDataPathFileName}	{V:BatchDataPathFileName}		
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:AnalystName}	{V:AnalystName}
Report Generation Time	{Expression...}	{C:ReportGeneratorName}	{Expression...}
{L:Calibration Last Update}	{V:CalibrationLastUpdatedTimeStamp}	{C:BatchState}	{V:BatchState}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}	{C:ReportQuantVersion}	{Expression...}
{V:CompoundName}	{C:RelativeStandardError}	{V:RelativeStandardError}	
CompoundCalibration			
{C:CalibrationSTDPPathName}	{C:CalibratiC:LevelName:LevelEnabl		
{V:CalibrationSTDPPathName}	{V:CalibrationV:LevelNameV:LevelEnable		
		{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

2 Report Template Workflows

Add a Calibration Results table

- 22 For the cells in the fifth column, select **Calibration** and **LevelConcentration** for the **Field Caption** of the cell in the first row, and **Calibration** and **LevelConcentration** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and set the horizontal alignment of both cells to **Right**. Click **OK**.

{L:Report Title}			
{C:BatchDataPathFileName}	{V:BatchDataPathFileName}		
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:AnalystName}	{V:AnalystName}
Report Generation Time	{Expression...}	{C:ReportGeneratorName}	{Expression...}
{L:Calibration Last Update}	{V:CalibrationLastUpdatedTimeStamp}	{C:BatchState}	{V:BatchState}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}	{C:ReportQuantVersion}	{Expression...}
{V:CompoundName}	{C:RelativeStandardError}	{V:RelativeStandardError}	
CompoundCalibration			
{C:CalibrationSTDPPathName}	{C:CalibratC:LevelName::LevelEnableConcentration}		
{V:CalibrationSTDPPathName}	{V:CalibrationV:LevelNameV:LevelEnableConcentration}		
		{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

- 23 To use the same format in the {V:LevelConcentration} cell as is found in the Quantitative Analysis software, double-click **Field Format** in the properties pane, and select **Calibration** from the **Binding Name** list, and select **LevelConcentration** from the **Field Name** list.

Textbox14

Field Format

Binding Name: Calibration

Field Name: LevelConcentration

Field Name

2 Report Template Workflows

Add a Calibration Results table

- 24 For the cells in the sixth column, select **Calibration** and **LevelResponse** for the **Field Caption** of the cell in the first row, and **Calibration** and **LevelResponse** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and set the horizontal alignment of both cells to **Right**.

The screenshot shows a report template with a table containing various fields. The properties pane on the right is open, showing the 'Field Format' settings for a selected cell. The 'Field Format' is set to 'Calibration-LevelResponse' and the 'Field Name' is also set to 'LevelResponse'. The 'Field Caption' is also set to 'Calibration-LevelResponse'.

- 25 Select the **{V:LevelResponse}** cell, double-click **Field Format** in the Properties pane, and select **Calibration** from the **Binding Name** list and **LevelResponse** from the **Field Name** list.

The screenshot shows the 'Field Format' section of the Properties pane. The 'Field Format' is set to 'Calibration-LevelResponse' and the 'Field Name' is also set to 'LevelResponse'. The 'Binding Name' is set to 'Calibration'.

2 Report Template Workflows

Add a Calibration Results table

- 26 For the cells in the seventh column, select **Calibration** and **LevelResponseFactor** for the **Field Caption** of the cell in the first row, and **Calibration** and **LevelResponseFactor** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and set the horizontal alignment of both cells to **Right**.

The screenshot shows a report template with a table containing various data points. The table has columns for different parameters and their values. The Properties pane on the right is open for a cell, showing the 'Field Caption' and 'Field Value' properties. The 'Field Caption' is set to 'Calibration-LevelResponseFactor' and the 'Field Value' is set to 'Textbox18'. The 'Field Format' is also set to 'Calibration-LevelResponseFactor'.

- 27 Select the **{V:LevelResponseFactor}** cell, double-click **Field Format** in the Properties pane, and select **Calibration** from the **Binding Name** list and **LevelResponseFactor** from the **Field Name** list.

The screenshot shows the Properties pane for a cell. The 'Field Format' section is expanded, and the 'Binding Name' is set to 'Calibration' and the 'Field Name' is set to 'LevelResponseFactor'. The 'General' section is also visible, showing the ID as 'Textbox18' and the Type as 'Text box'.

2 Report Template Workflows

Add a Calibration Results table

- 28 For the cells in the eighth column, select **Calibration** and **LevelRSD** for the **Field Caption** of the cell in the first row, and **Calibration** and **LevelRSD** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and set the horizontal alignment of both cells to **Right**.

The screenshot shows a report template with a table containing various fields. The table has columns for different parameters and their values. The Properties pane on the right is open for a text box, showing settings for Borders, Padding, Content, Field Caption, Field Value, Text, Fill, Background Color, Font, and Format. The Format field is set to F1.

- 29 Select the **{V:LevelRSD}** cell, and enter **F1** in the **Format** field in the Properties pane.

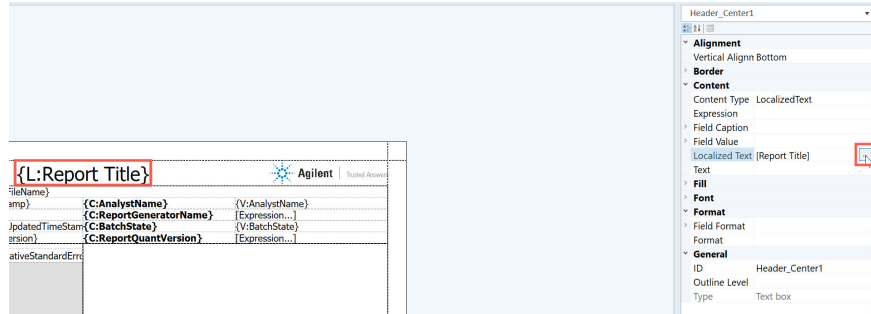
The screenshot shows the Properties pane for a text box. The Format field is highlighted with a red box and contains the text 'F1'. Other fields like Alignment, Border, Content, Fill, Font, and General are also visible.

2 Report Template Workflows

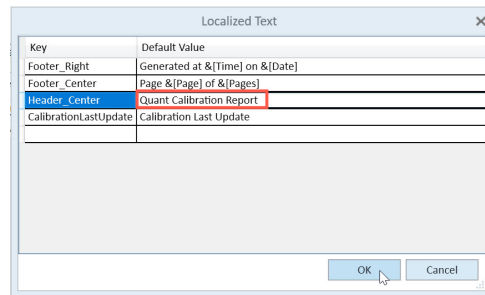
Add a report title

Add a report title

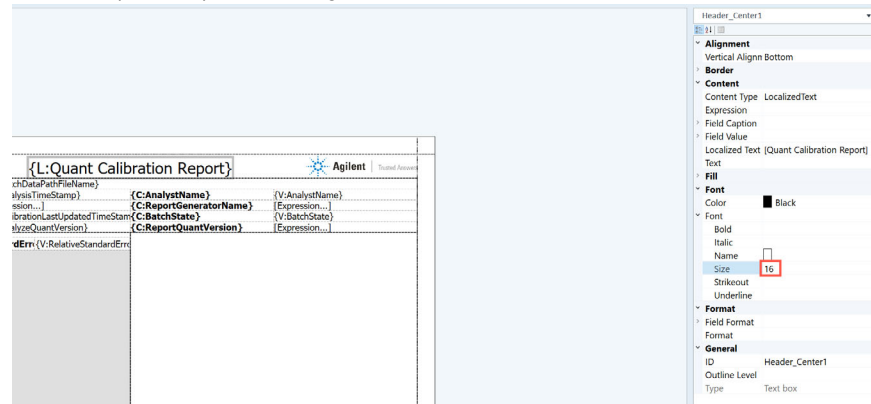
- 1 Select the **{L:Report Title}** cell in the page header. Click **Localized Text** in the Properties pane, then click the ellipsis.



- 2 Enter **Quant Calibration Report** in the **Default Value** column for **Header Center**. Click **OK**.



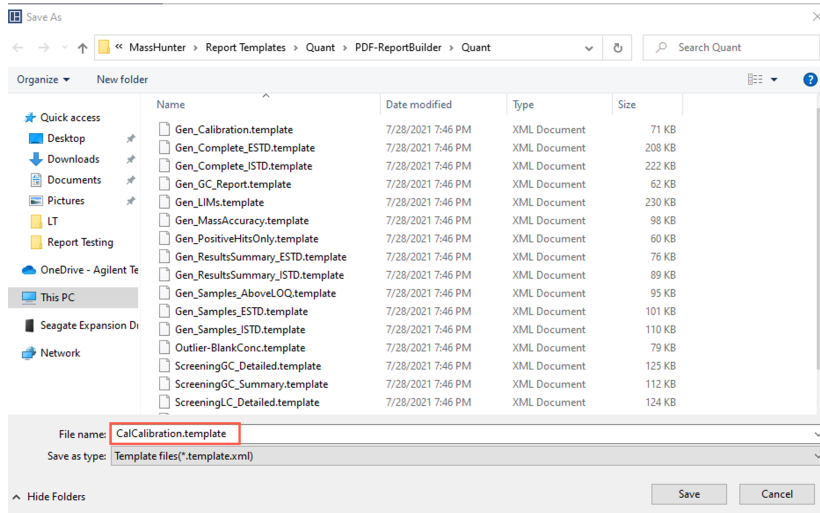
- 3 In the Properties pane, change the font size to **16**.



2 Report Template Workflows

Add a report title

- 4 Click **File > Save As**, and save the template as **CalCalibration.template.xml**.

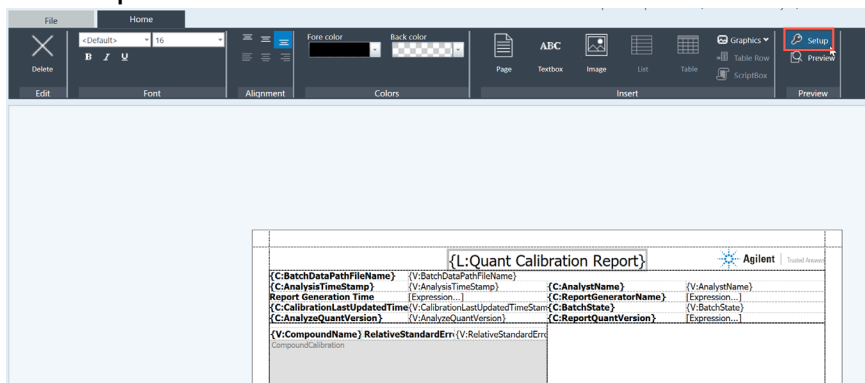


2 Report Template Workflows

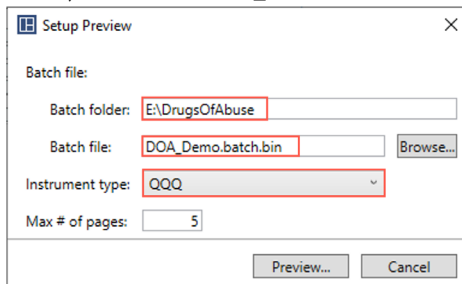
Preview the report template

Preview the report template

- 1 Click **Setup**.



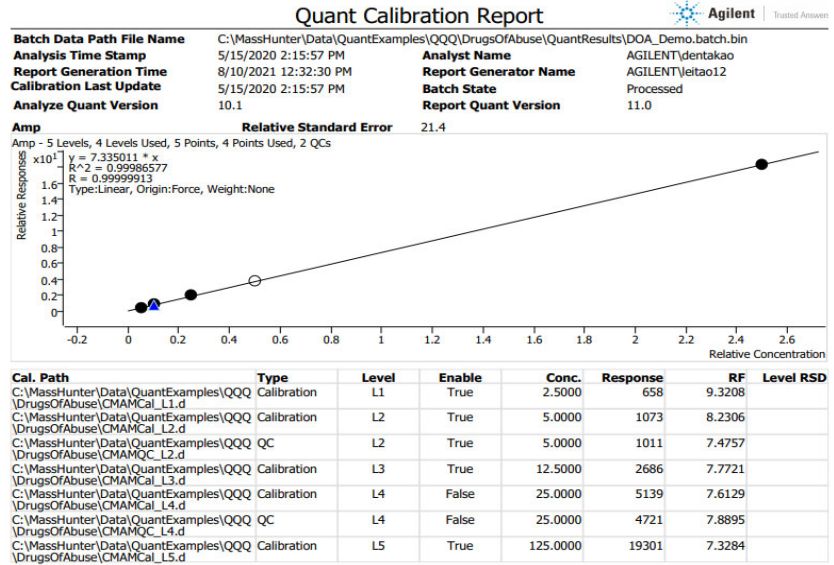
- 2 Click **browse** in the Setup Preview window, navigate to the **DrugsOfAbuse** folder, and select **DOA_Demo.batch.bin**. Select **QQQ** for **Instrument type**.



2 Report Template Workflows

Preview the report template

3 Click **Preview**. A PDF report is generated using the selected batch file.



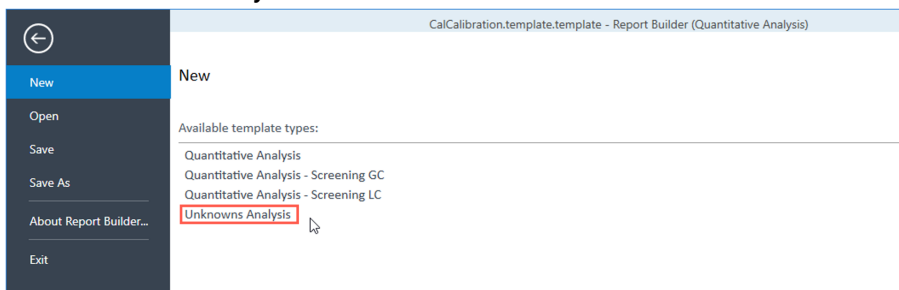
Unknowns Analysis - Area Percent Report Template

For this example, you will create a Calibration Report by completing the following procedures.

- 1 **Create a new Unknowns Analysis report template** on page 50
- 2 **Insert a list** on page 51
- 3 **Add a Sample Information table with a graph** on page 53
- 4 **Add a Compound Results table** on page 61
- 5 **Add a report title** on page 91
- 6 **Generate a report** on page 93

Create a new Unknowns Analysis report template

To create a new Unknowns Analysis report template, select **File > New**, and select **Unknowns Analysis**.

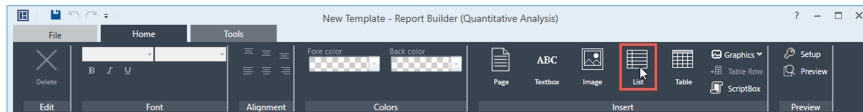


2 Report Template Workflows

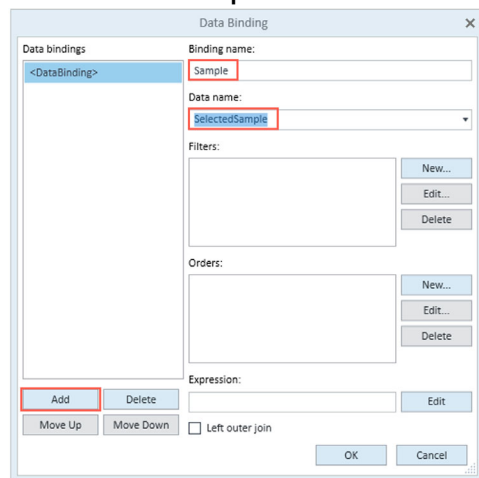
Insert a list

Insert a list

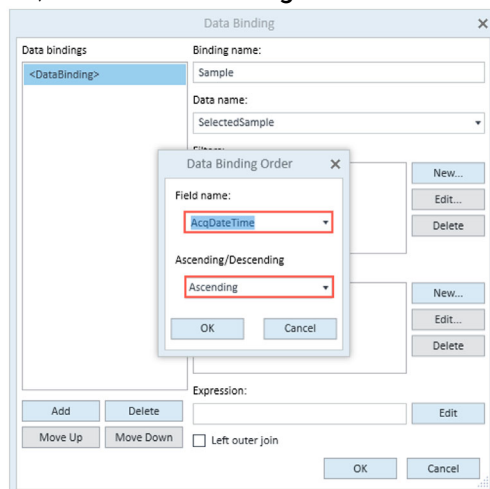
1 Click **List**.



2 Under **Data bindings**, click **Add**. Enter **Sample** in the **Binding name** field, and select **SelectedSample** from the **Data name** list.



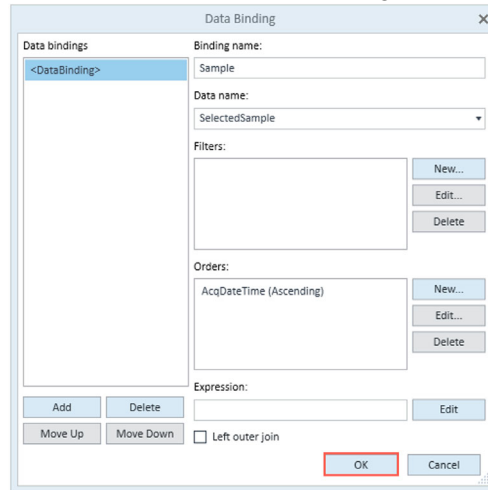
3 In the **Orders** section, click **New**. Select **AcqDateTime** from the **Field name** list, and select **Ascending** under the **Ascending/Descending** list. Click **OK**.



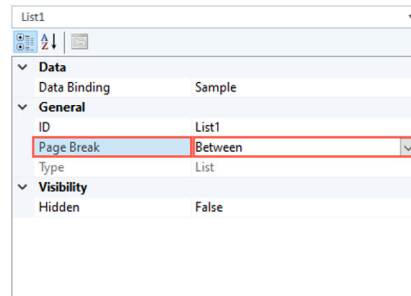
2 Report Template Workflows

Insert a list

- 4 Click **OK** to close the **Data Binding** window.



- 5 Select **Page Break** in the Properties pane, and select **Between** from the list.

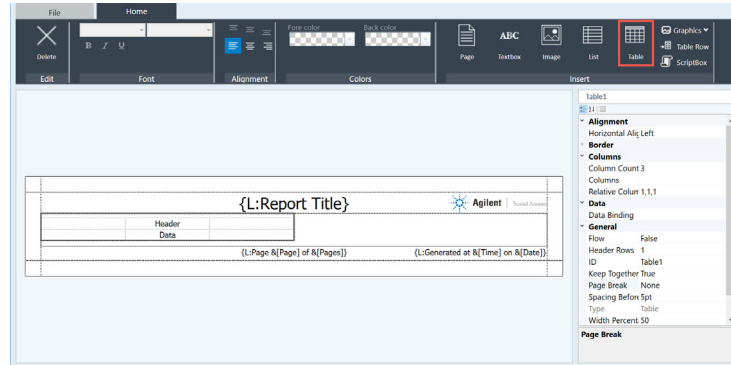


2 Report Template Workflows

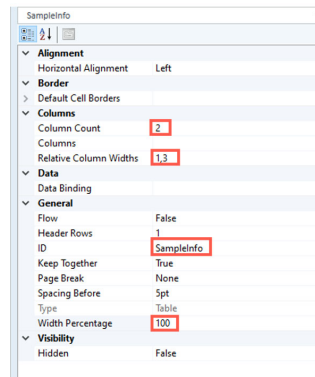
Add a Sample Information table with a graph

Add a Sample Information table with a graph

- 1 Select **List1** from the drop-down menu at the top of the Properties pane. Click **Table**. A new table (Table1) appears in the report.



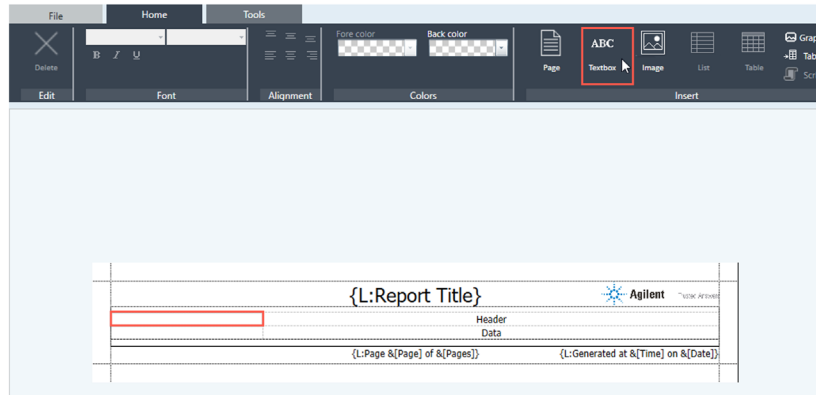
- 2 In the Properties pane, make the following changes:
 - Change the **Column Count** to 2
 - Change the **Relative Column Widths** to 1,3 to resize the columns in relation to the table width
 - Enter **SampleInfo** in the **ID** field to update the table name
 - Enter 100 in the **Width Percentage** field to resize the table width.



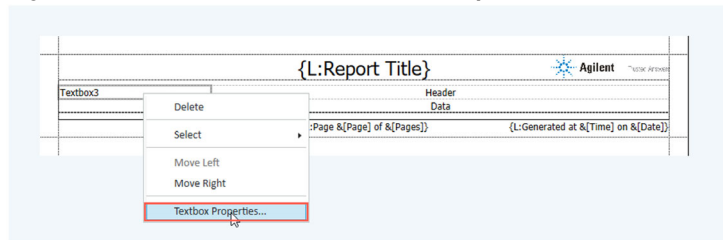
2 Report Template Workflows

Add a Sample Information table with a graph

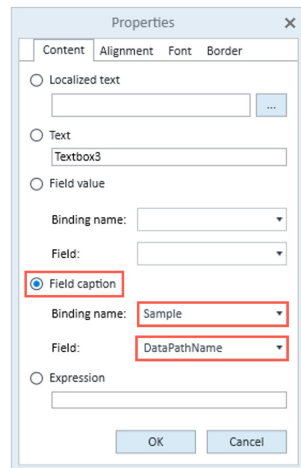
- 3 To edit a table cell, select the empty cell to the left of the **Header** cell, and click **Textbox**.



- 4 Right-click the cell and select **Textbox Properties**.



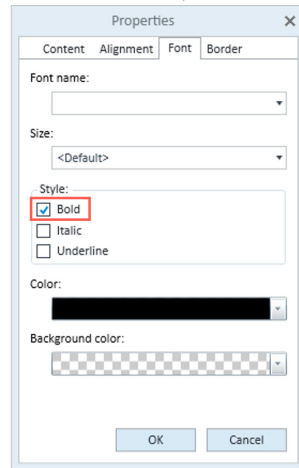
- 5 Select **Field caption**, then select **Sample** from the **Binding name** list, and select **DataPathName** from the **Field** list.



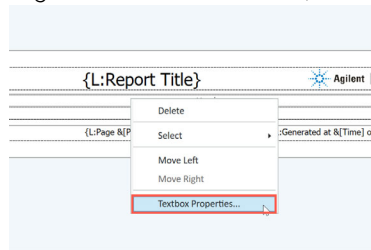
2 Report Template Workflows

Add a Sample Information table with a graph

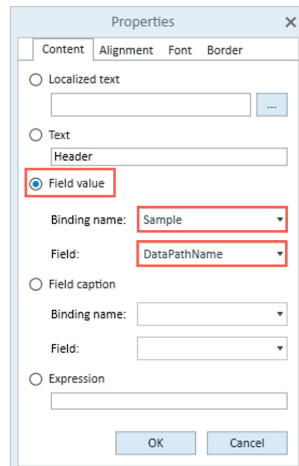
- 6 On the **Font** tab, select **Bold**. Click **OK**.



- 7 Right-click the **Header** cell, and select **Textbox Properties**.



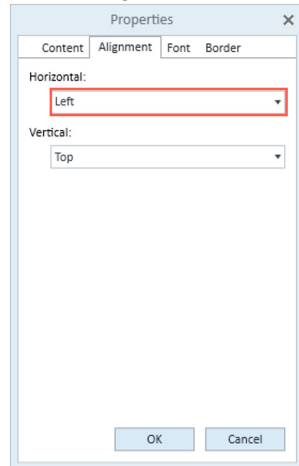
- 8 Select **Field Value**, select **Sample** from the **Binding name** list, and **DataPathName** from the **Field** list.



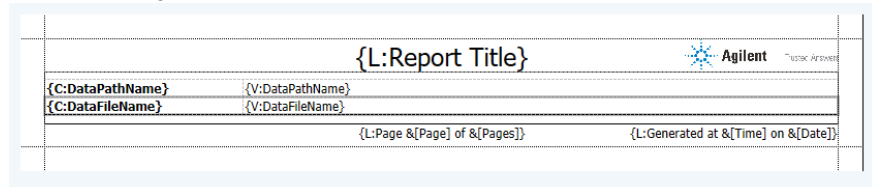
2 Report Template Workflows

Add a Sample Information table with a graph

- 9 On the **Alignment** tab, select **Left** from the **Horizontal** list. Click **OK**.



- 10 For the cells in the second row, use the instructions in the previous steps to select **Sample** and **DataFileName** for the **Field Caption** of the cell in the first column, and **Sample** and **DataFileName** for the **Field Value** of the cell in the second column. Bold the text of the cell in the first column, and set the horizontal alignment of both cells to **Left**. Click **OK**.



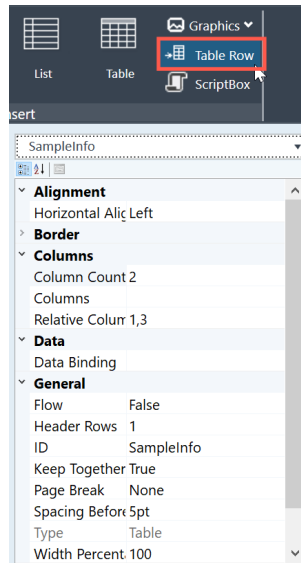
The image shows a report template layout. At the top right is the Agilent logo and 'Agilent' text. Below it is a table with the following content:

{L:Report Title}	
{C:DataPathName}	{V:DataPathName}
{C:DataFileName}	{V:DataFileName}
{L:Page &[Page] of &[Pages]}	
{L:Generated at &[Time] on &[Date]}	

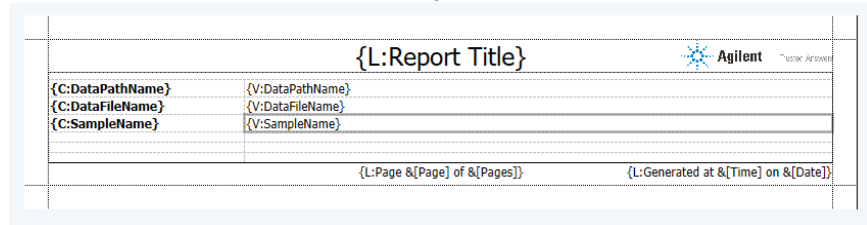
2 Report Template Workflows

Add a Sample Information table with a graph

- 11 To add rows to the table, select **SampleInfo** from the list at the top of the Properties pane, and click **Table Row**. Click **Table Row** 3 more times to add a total of 4 rows.



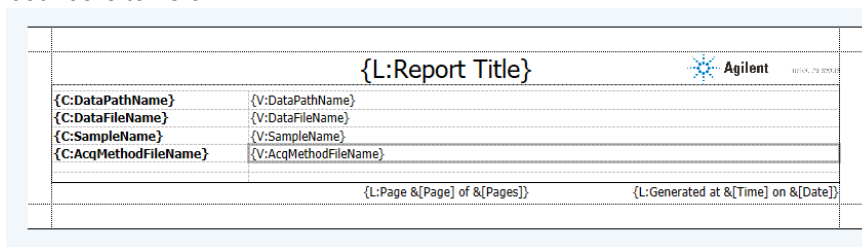
- 12 For the cells in the third row, select **Sample** and **SampleName** for the **Field Caption** of the cell in the first column, and **Sample** and **SampleName** for the **Field Value** of the cell in the second column. Bold the text of the cell in the first column, and set the horizontal alignment of both cells to **Left**.



2 Report Template Workflows

Add a Sample Information table with a graph

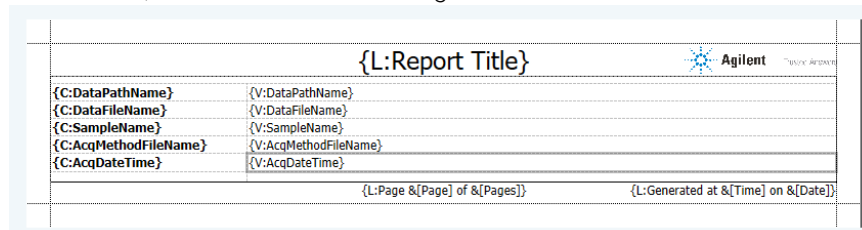
- 13 For the cells in the fourth row, select **Sample** and **AcqMethodFileName** for the **Field Caption** of the cell in the first column, and **Sample** and **AcqMethodFileName** for the **Field Value** of the cell in the second column. Bold the text of the cell in the first column, and set the horizontal alignment of both cells to **Left**.



The screenshot shows a report template with a table containing the following data:

{L:Report Title}		Agilent	10/14, 2015
{C:DataPathName}	{V:DataPathName}		
{C:DataFileName}	{V:DataFileName}		
{C:SampleName}	{V:SampleName}		
{C:AcqMethodFileName}	{V:AcqMethodFileName}		
		{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

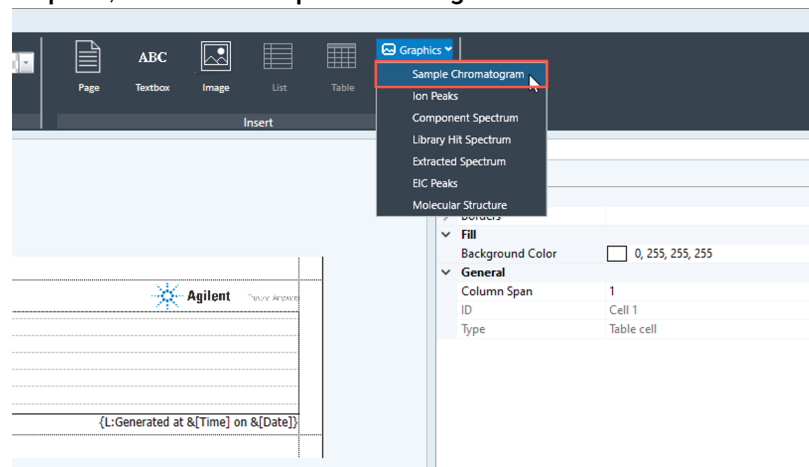
- 14 For the cells in the fifth row, select **Sample** and **AcqDateTime** for the **Field Caption** of the cell in the first column, and **Sample** and **AcqDateTime** for the **Field Value** of the cell in the second column. Bold the text of the cell in the first column, and set the horizontal alignment of both cells to **Left**.



The screenshot shows a report template with a table containing the following data:

{L:Report Title}		Agilent	10/14, 2015
{C:DataPathName}	{V:DataPathName}		
{C:DataFileName}	{V:DataFileName}		
{C:SampleName}	{V:SampleName}		
{C:AcqMethodFileName}	{V:AcqMethodFileName}		
{C:AcqDateTime}	{V:AcqDateTime}		
		{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

- 15 To add a graphic, select the empty cell below **{C:AcqDateTime}**, click **Graphics**, and select **Sample Chromatogram**.



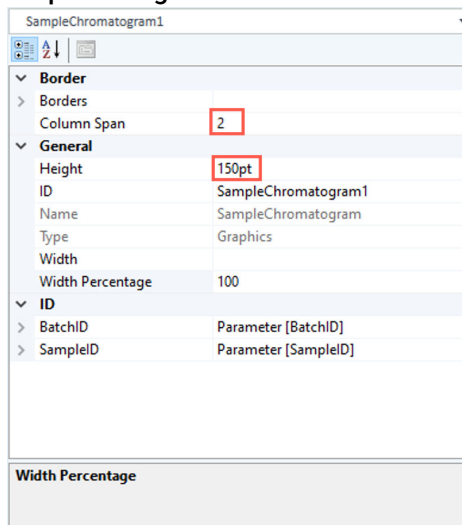
The screenshot shows the Report Builder interface. The **Graphics** menu is open, and **Sample Chromatogram** is selected. The table below the menu contains the following data:

{L:Report Title}		Agilent	10/14, 2015
{C:DataPathName}	{V:DataPathName}		
{C:DataFileName}	{V:DataFileName}		
{C:SampleName}	{V:SampleName}		
{C:AcqMethodFileName}	{V:AcqMethodFileName}		
{C:AcqDateTime}	{V:AcqDateTime}		
		{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

2 Report Template Workflows

Add a Sample Information table with a graph

- 16 To resize the graphic, in the Properties pane, enter **2** for **Column Span** and **150pt** for **Height**.



- 17 ID fields like BatchID and SampleID have system-assigned numbers which help to rapidly locate and link to unique table rows in a Quantitative Analysis dataset. To edit the ID fields, expand the ID section in the Properties pane. Enter the values for each ID:

- BatchID
 - **Field Value:** Sample-BatchID
 - **Binding Name:** Sample
 - **Field Name:** BatchID
- SampleID
 - **Field Value:** Sample-SampleID
 - **Binding Name:** Sample

2 Report Template Workflows

Add a Sample Information table with a graph

- **Field Name:** SampleID

The screenshot shows a configuration window for a report element named 'SampleChromatogram1'. The window is divided into two main sections: a top section for general properties and a bottom section for field configuration.

General Properties:

Height	150pt
ID	SampleChromatogram1
Name	SampleChromatogram
Type	Graphics
Width	
Width Percentage	100

Field Configuration:

- Expanded **ID** folder:
 - Expanded **BatchID** folder:
 - Field Value: Parameter [BatchID]
 - Field Value: Sample-BatchID
 - Binding Name: Sample
 - Field Name: BatchID
 - Value:
 - Expanded **SampleID** folder:
 - Field Value: Parameter [SampleID]
 - Field Value: Sample-SampleID
 - Binding Name: Sample
 - Field Name: SampleID
 - Value:

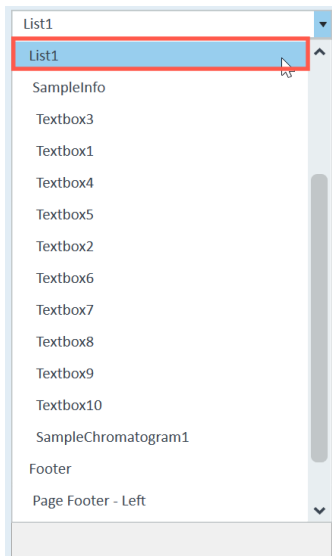
The 'Field Name' section at the bottom of the window is currently empty.

2 Report Template Workflows

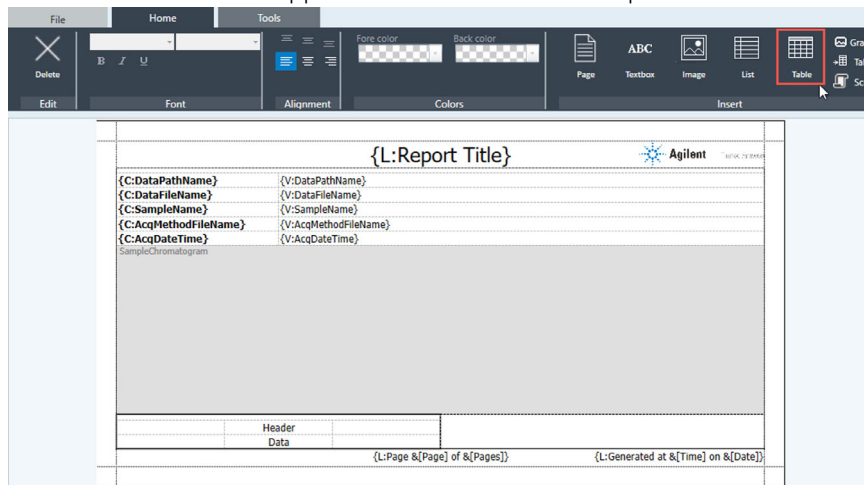
Add a Compound Results table

Add a Compound Results table

- 1 Select **List1** from the list at the top of the Properties pane.



- 2 Click **Table**. A new table appears at the bottom of the report.

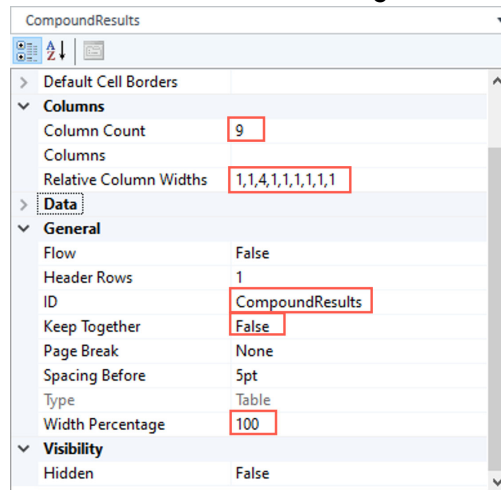


- 3 In the properties pane,
 - change the **Column Count** to 9
 - change the **Relative Column Width** to 1,1,4,1,1,1,1,1,1 to resize the columns in relation to the table width

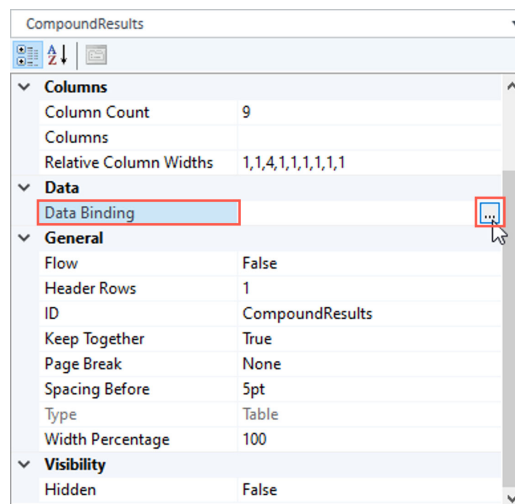
2 Report Template Workflows

Add a Compound Results table

- enter CompoundResults in the **ID** field to update the table name
- select False for **Keep Together**
- enter 100 in the **Width Percentage** field to resize the table width



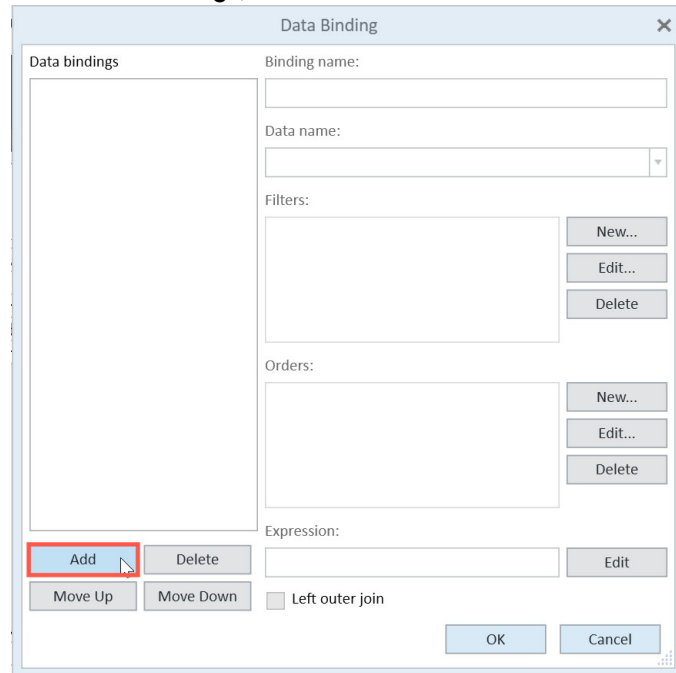
- 4 To add binding data to the table, click **Data Binding** in the Properties pane, then click the ellipsis.



2 Report Template Workflows

Add a Compound Results table

5 Under **Data Bindings**, click **Add**.



The screenshot shows a dialog box titled "Data Binding" with a close button (X) in the top right corner. The dialog is divided into several sections:

- Data bindings:** A large empty list box on the left side.
- Binding name:** A text input field.
- Data name:** A dropdown menu.
- Filters:** A list box with three buttons to its right: "New...", "Edit...", and "Delete".
- Orders:** A list box with three buttons to its right: "New...", "Edit...", and "Delete".
- Expression:** A text input field with an "Edit" button to its right.
- Left outer join:** A checkbox.
- Bottom controls:** A row of buttons: "Add" (highlighted with a red box and a mouse cursor), "Delete", "Move Up", "Move Down", "OK", and "Cancel".

2 Report Template Workflows

Add a Compound Results table

- 6 Enter **Component** for the **Binding name** and select **Component** from the **Data name** list.

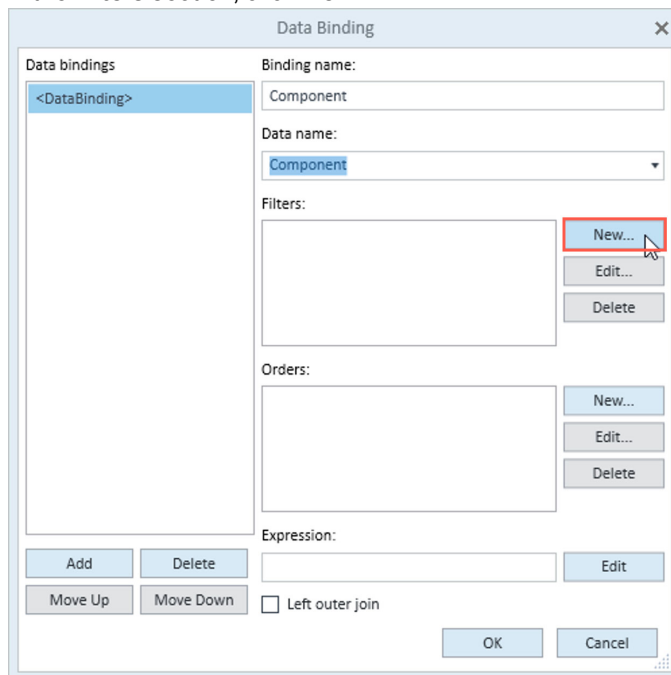
The screenshot shows the 'Data Binding' dialog box with the following fields and controls:

- Data bindings:** A list containing '<DataBinding>'. Below the list are buttons for 'Add', 'Delete', 'Move Up', and 'Move Down'.
- Binding name:** A text field containing 'Component'.
- Data name:** A dropdown menu with 'Component' selected.
- Filters:** An empty list box with buttons for 'New...', 'Edit...', and 'Delete'.
- Orders:** An empty list box with buttons for 'New...', 'Edit...', and 'Delete'.
- Expression:** An empty text field with an 'Edit' button.
- Left outer join:** A checkbox that is currently unchecked.
- Buttons:** 'OK' and 'Cancel' buttons at the bottom right.

2 Report Template Workflows

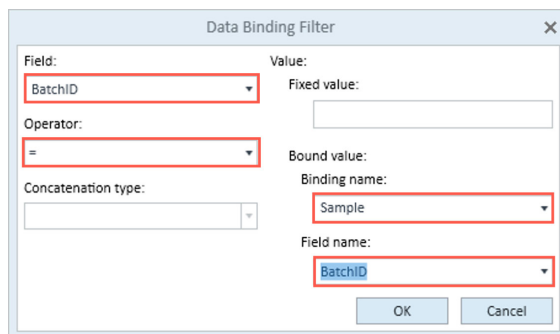
Add a Compound Results table

- 7 In the **Filters** section, click **New**.



The screenshot shows the 'Data Binding' dialog box. On the left, there is a list of data bindings with '<DataBinding>' selected. On the right, there are fields for 'Binding name:' (Component), 'Data name:' (Component), and 'Filters:'. The 'Filters:' section has a red box around the 'New...' button. Below it are 'Orders:' and 'Expression:' sections. At the bottom, there are 'Add', 'Delete', 'Move Up', 'Move Down', 'Left outer join', 'OK', and 'Cancel' buttons.

- 8 Select **BatchID** from the **Field** list, and **=** from the **Operator** list. Select **Sample** from the **Binding name** list, and select **BatchID** from the **Field name** list. Click **OK**.



The screenshot shows the 'Data Binding Filter' dialog box. It has fields for 'Field:' (BatchID), 'Operator:' (=), 'Concatenation type:', 'Value:', 'Fixed value:', 'Bound value:', 'Binding name:' (Sample), and 'Field name:' (BatchID). The 'Field', 'Operator', 'Binding name', and 'Field name' dropdowns are highlighted with red boxes. At the bottom, there are 'OK' and 'Cancel' buttons.

2 Report Template Workflows

Add a Compound Results table

9 In the **filters** section, click **New**.

The screenshot shows the 'Data Binding' dialog box. On the left, there is a list of data bindings with '<DataBinding>' selected. The right side contains configuration fields: 'Binding name' (Component), 'Data name' (Component), and 'Filters' (<BatchID> = <Sample:BatchID>). The 'New...' button in the 'Filters' section is highlighted with a red box. Below the filters are 'Orders' and 'Expression' sections, each with 'New...', 'Edit...', and 'Delete' buttons. At the bottom, there are 'Add', 'Delete', 'Move Up', 'Move Down', and 'Left outer join' options, along with 'OK' and 'Cancel' buttons.

10 Select **SampleID** from the **Field** list, **Sample** from the **Binding Name** list, and **SampleID** from the **Field Name** list.

The screenshot shows the 'Data Binding Filter' dialog box. It has several fields: 'Field' (SampleID), 'Value' (empty), 'Operator' (=), 'Concatenation type' (AND), 'Fixed value' (empty), 'Bound value' (empty), 'Binding name' (Sample), and 'Field name' (SampleID). The 'Field', 'Binding name', and 'Field name' dropdowns are highlighted with red boxes. 'OK' and 'Cancel' buttons are at the bottom.

2 Report Template Workflows

Add a Compound Results table

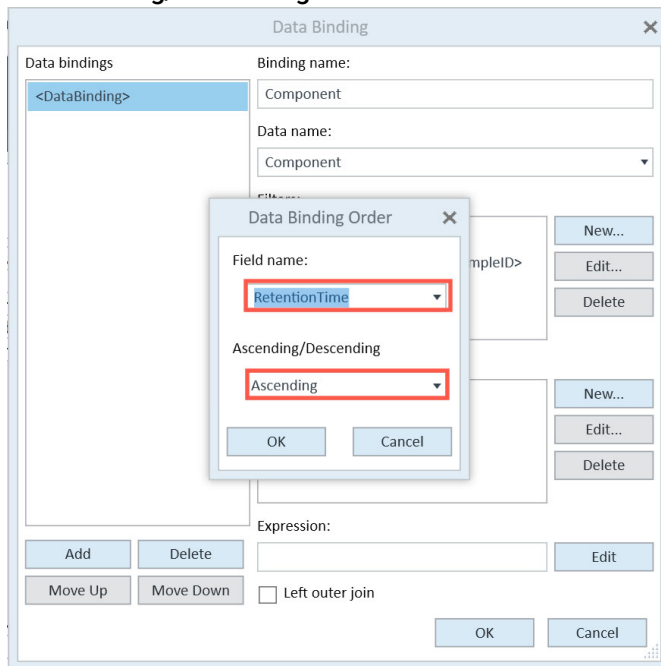
11 In the **Orders** section, click **New**.

The screenshot shows the 'Data Binding' dialog box. The 'Data bindings' list on the left contains a single entry '<DataBinding>'. The 'Binding name:' field is set to 'Component'. The 'Data name:' dropdown is also set to 'Component'. The 'Filters:' section contains two filter expressions: '<BatchID> = <Sample:BatchID>' and 'AND <SampleID> = <Sample:SampleID>'. The 'Orders:' section is currently empty, and the 'New...' button next to it is highlighted with a red rectangular box. At the bottom of the dialog, there are buttons for 'Add', 'Delete', 'Move Up', 'Move Down', 'Expression:', 'Left outer join' (checkbox), 'Edit', 'OK', and 'Cancel'.

2 Report Template Workflows

Add a Compound Results table

- 12 Select **RetentionTime** from the **Field name** list, and select **Ascending** from the **Ascending/Descending** list. Click **OK**.



2 Report Template Workflows

Add a Compound Results table

13 Under **Data Bindings**, click **Add**.

The screenshot shows the 'Data Binding' dialog box. The 'Data bindings' list on the left contains a single entry '<DataBinding>'. The 'Binding name:' field is set to 'Component'. The 'Data name:' dropdown is also set to 'Component'. The 'Filters:' section contains two filter expressions: '<BatchID> = <Sample:BatchID>' and 'AND <SampleID> = <Sample:SampleID>'. The 'Orders:' section contains one order expression: 'RetentionTime (Ascending)'. At the bottom, the 'Add' button is highlighted with a red box and a mouse cursor. Other buttons include 'Delete', 'Move Up', 'Move Down', 'Expression:', 'Left outer join' (checkbox), 'OK', and 'Cancel'.

2 Report Template Workflows

Add a Compound Results table

14 Enter **Hit** in the **Binding name** field, and select **Hit** from the **Data name** list.

The screenshot shows the 'Data Binding' dialog box. On the left, a list of 'Data bindings' includes 'Component' and '<DataBinding>'. The 'Binding name' field is set to 'Hit'. The 'Data name' dropdown menu is open, showing 'Hit' selected. Below this are sections for 'Filters' and 'Orders', each with 'New...', 'Edit...', and 'Delete' buttons. At the bottom, there are buttons for 'Add', 'Delete', 'Move Up', 'Move Down', 'Expression', and 'Left outer join', along with 'OK' and 'Cancel' buttons.

15 Add a data binding filter for each of the following configurations:

- a Select **BatchID** from the **Field** list, select **=** from the **Operator** list, select **Sample** from the **Binding name** list, and select **BatchID** from the **Field name** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'BatchID', the 'Operator' dropdown is set to '=', the 'Binding name' dropdown is set to 'Sample', and the 'Field name' dropdown is set to 'BatchID'. There are also fields for 'Value', 'Fixed value', 'Bound value', and 'Concatenation type'. At the bottom, there are 'OK' and 'Cancel' buttons.

2 Report Template Workflows

Add a Compound Results table

- b Select **SampleID** from the **Field** list, select **=** from the **Operator** list, select **AND** from the **Concatenation type** list, select **Sample** from the **Binding name** list, and select **SampleID** from the **Field name** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'SampleID', the 'Operator' dropdown is set to '=', the 'Concatenation type' dropdown is set to 'AND', the 'Binding name' dropdown is set to 'Sample', and the 'Field name' dropdown is set to 'SampleID'. The 'Fixed value' and 'Bound value' fields are empty. The 'OK' and 'Cancel' buttons are at the bottom right.

- c Select **ComponentID** from the **Field** list, select **=** from the **Operator** list, select **AND** from the **Concatenation type** list, select **Component** from the **Binding name** list, and select **ComponentID** from the **Field name** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'ComponentID', the 'Operator' dropdown is set to '=', the 'Concatenation type' dropdown is set to 'AND', the 'Binding name' dropdown is set to 'Component', and the 'Field name' dropdown is set to 'ComponentID'. The 'Fixed value' and 'Bound value' fields are empty. The 'OK' and 'Cancel' buttons are at the bottom right.

2 Report Template Workflows

Add a Compound Results table

- d Select **HitID** from the **Field** list, select **=** from the **Operator** list, select **AND** from the **Concatenation type** list, select **Component** from the **Binding name** list, and select **PrimaryHitID** from the **Field name** list.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'HitID', the 'Operator' dropdown is set to '=', the 'Concatenation type' dropdown is set to 'AND', the 'Binding name' dropdown is set to 'Component', and the 'Field name' dropdown is set to 'PrimaryHitID'. The 'Fixed value' field is empty. The 'OK' and 'Cancel' buttons are at the bottom right.

- e Select **Visible** from the **Field** list, select **=** from the **Operator** list, select **AND** from the **Concatenation type** list, and enter **True** in the **Fixed value** field.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' dropdown is set to 'Visible', the 'Operator' dropdown is set to '=', the 'Concatenation type' dropdown is set to 'AND', and the 'Fixed value' field contains 'True'. The 'Binding name' and 'Field name' fields are empty. The 'OK' and 'Cancel' buttons are at the bottom right.

2 Report Template Workflows

Add a Compound Results table

16 Under **Data Bindings**, select **Component**, and select **Left outer join**.

The screenshot shows the 'Data Binding' dialog box. On the left, under 'Data bindings', 'Component' is selected and highlighted with a red box. The right side of the dialog contains the following fields and controls:

- Binding name:** Component
- Data name:** Component
- Filters:** <BatchID> = <Sample:BatchID> AND <SampleID> = <Sample:SampleID>. Includes 'New...', 'Edit...', and 'Delete' buttons.
- Orders:** RetentionTime (Ascending). Includes 'New...', 'Edit...', and 'Delete' buttons.
- Expression:** (empty field) with 'Add', 'Delete', and 'Edit' buttons.
- Join type:** A checkbox labeled 'Left outer join' is checked and highlighted with a red box.
- Navigation:** 'Add', 'Delete', 'Move Up', and 'Move Down' buttons.
- Final Buttons:** 'OK' and 'Cancel' buttons.

2 Report Template Workflows

Add a Compound Results table

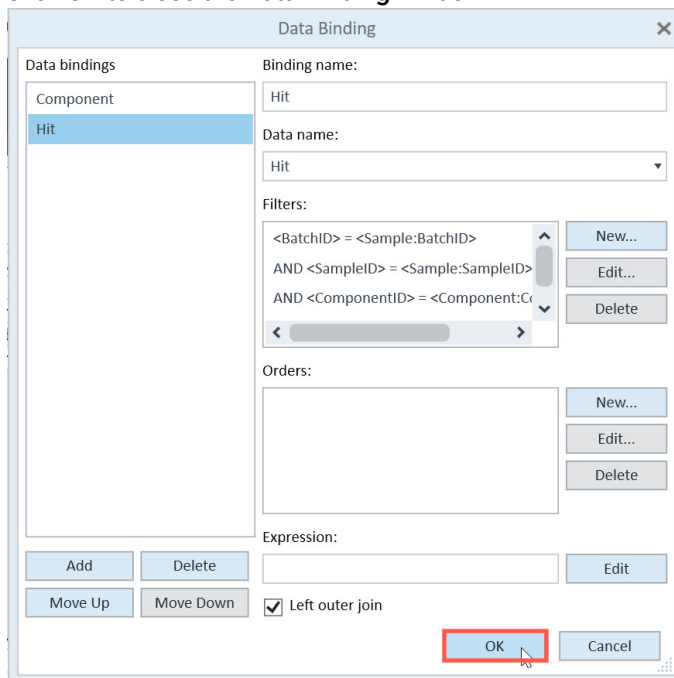
- 17 Under **Data Bindings**, select **Hit**, and select **Left outer join**. All components and matching hits are now included in the result table.

The screenshot shows the 'Data Binding' dialog box. On the left, under 'Data bindings', there is a table with two rows: 'Component' and 'Hit'. The 'Hit' row is highlighted with a red border. To the right of this table are fields for 'Binding name:' (containing 'Hit') and 'Data name:' (containing 'Hit'). Below these are 'Filters:' and 'Orders:' sections, each with a list of conditions and associated 'New...', 'Edit...', and 'Delete' buttons. At the bottom, there is an 'Expression:' field and a 'Left outer join' checkbox, which is checked and highlighted with a red border. Other buttons at the bottom include 'Add', 'Delete', 'Move Up', 'Move Down', 'OK', and 'Cancel'.

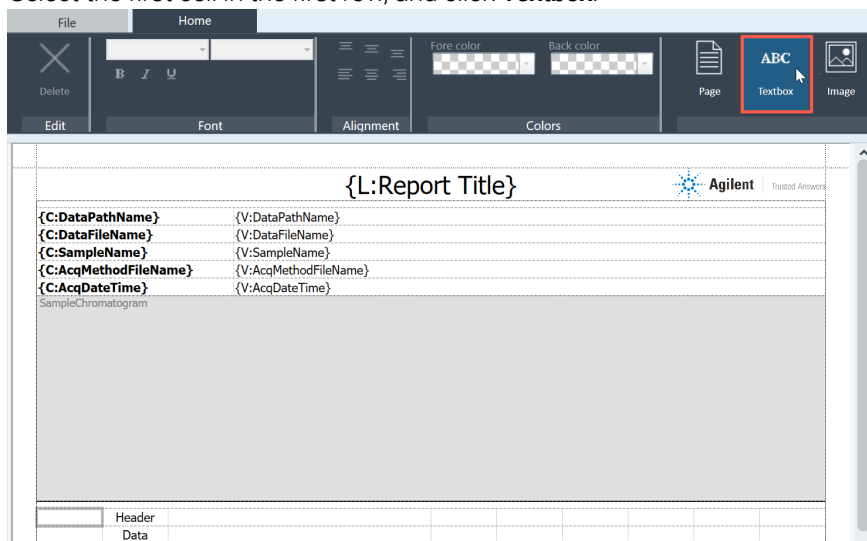
2 Report Template Workflows

Add a Compound Results table

18 Click **OK** to close the **Data Binding** window.



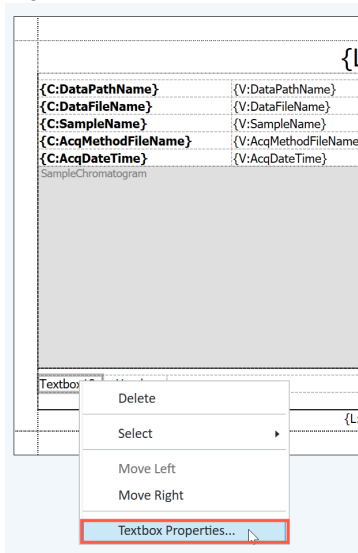
19 Select the first cell in the first row, and click **Textbox**.



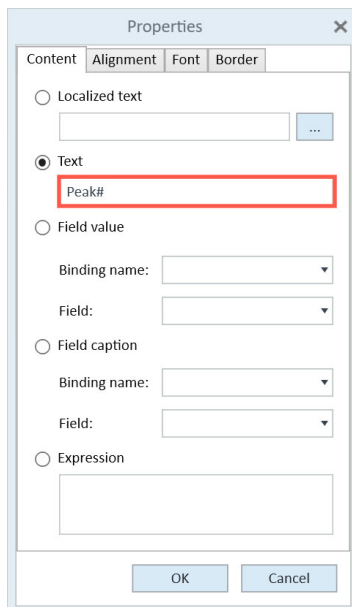
2 Report Template Workflows

Add a Compound Results table

20 Right-click the cell, and select **Textbox Properties**.



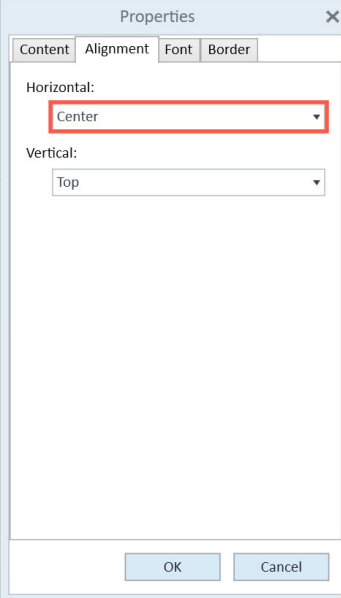
21 Enter **Peak#** in the **Text** field.



2 Report Template Workflows

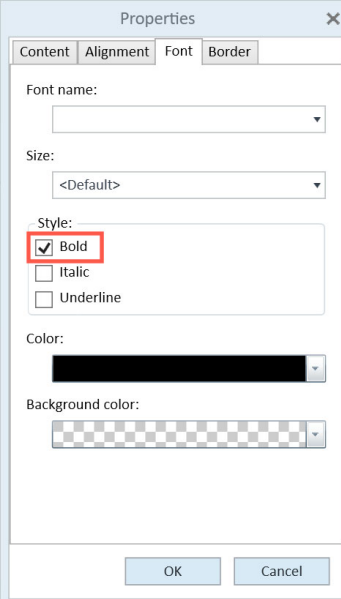
Add a Compound Results table

22 On the **Alignment** tab, select **Center** from the **Horizontal** list.



The screenshot shows the 'Properties' dialog box with the 'Alignment' tab selected. The 'Horizontal' dropdown menu is set to 'Center' and is highlighted with a red rectangle. The 'Vertical' dropdown menu is set to 'Top'. The 'Content', 'Font', and 'Border' tabs are also visible at the top of the dialog. 'OK' and 'Cancel' buttons are at the bottom.

23 On the **Font** tab, select **Bold**. Click **OK**.

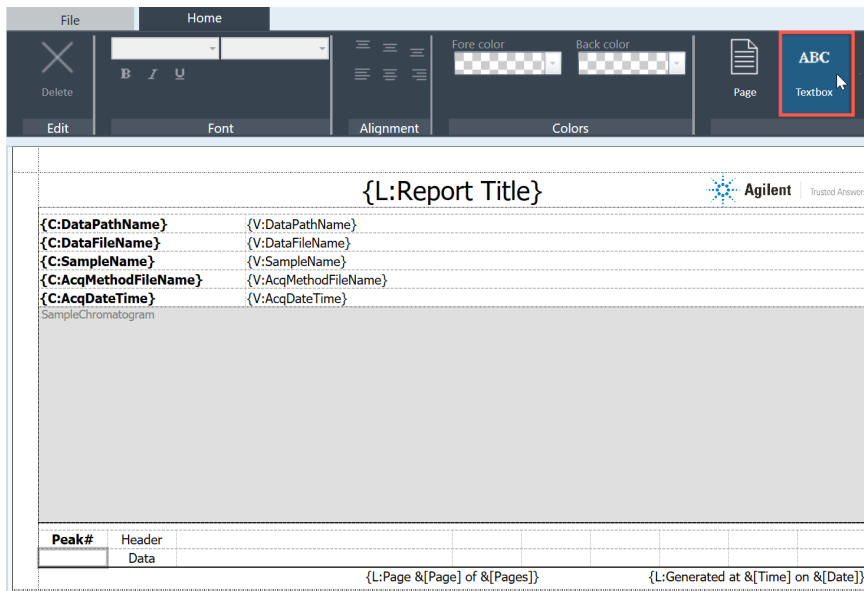


The screenshot shows the 'Properties' dialog box with the 'Font' tab selected. The 'Style' section has the 'Bold' checkbox checked and highlighted with a red rectangle. Other options like 'Italic' and 'Underline' are unchecked. The 'Font name' dropdown is empty, 'Size' is '<Default>', 'Color' is black, and 'Background color' is a checkerboard pattern. 'OK' and 'Cancel' buttons are at the bottom.

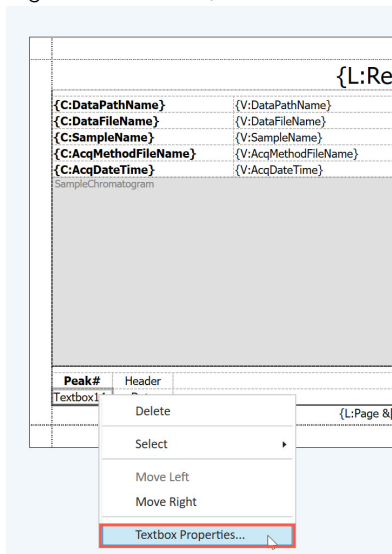
2 Report Template Workflows

Add a Compound Results table

24 Select the first cell in the second row and click **Textbox**.



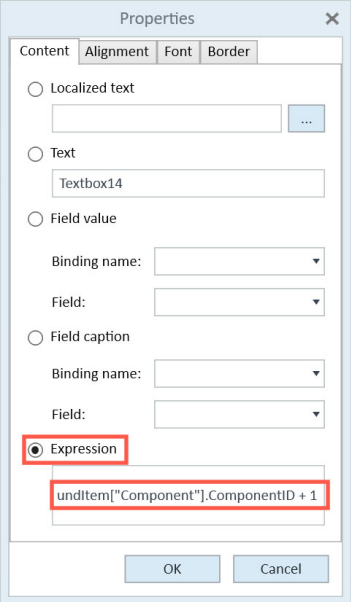
25 Right-click the cell, and select **Textbox Properties**.



2 Report Template Workflows

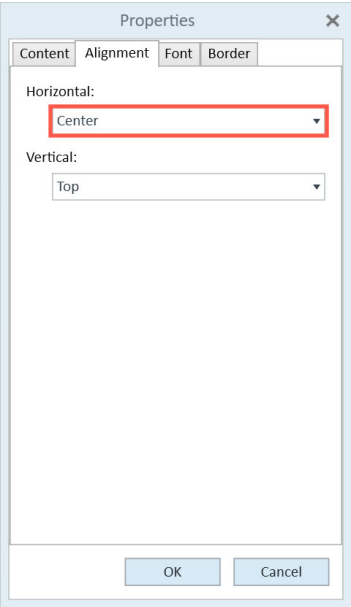
Add a Compound Results table

26 Select **Expression**, and enter **BoundItems["Component"].ComponentID + 1**.



The screenshot shows the 'Properties' dialog box with the 'Content' tab selected. The 'Expression' radio button is selected and highlighted with a red box. Below it, the text box contains the expression 'BoundItems["Component"].ComponentID + 1', which is also highlighted with a red box. The 'OK' and 'Cancel' buttons are visible at the bottom.

27 On the Alignment tab, select **Center** from the **Horizontal** list.

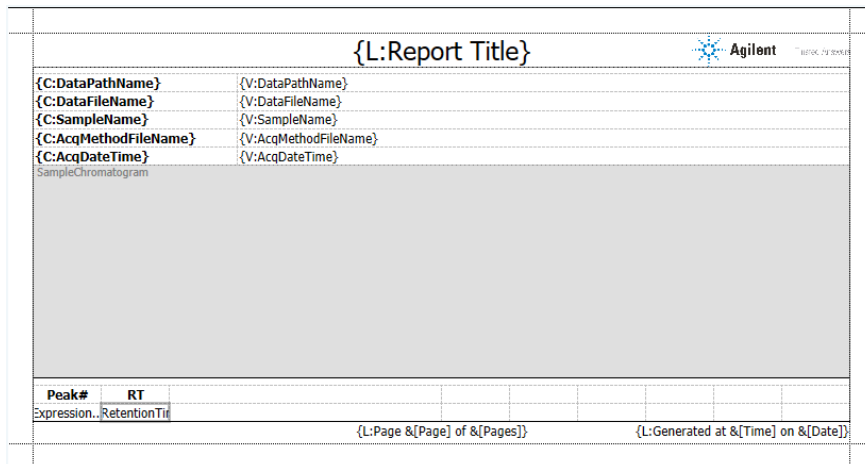


The screenshot shows the 'Properties' dialog box with the 'Alignment' tab selected. The 'Horizontal' dropdown menu is set to 'Center' and is highlighted with a red box. The 'Vertical' dropdown menu is set to 'Top'. The 'OK' and 'Cancel' buttons are visible at the bottom.

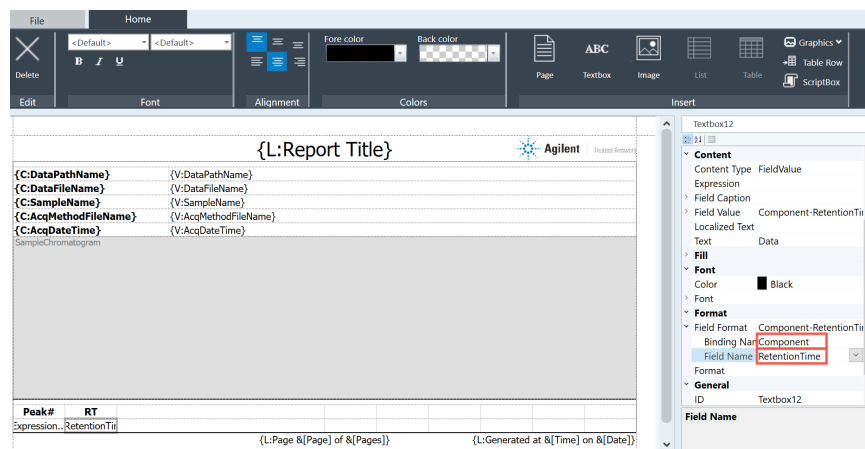
2 Report Template Workflows

Add a Compound Results table

- 28 For the cells in the second column, enter **RT** in the **Text** field for the cell in the first row and set the font to **Bold**. For the cell in the second row, select **Field value**, select **Component** from the **Binding name** list, and select **RetentionTime** from the **Field** list. Set the horizontal alignment of both cells to **Center**.



- 29 To use the same format in the **{V:RetentionTime}** cell as is found in the Quantitative Analysis software, double-click **Field Format** in the Properties pane, and select **Component** from the **Binding Name** list, and **RetentionTime** from the **Field Name** list.



2 Report Template Workflows

Add a Compound Results table

- 30 For the cells in the third column, use the instructions in the previous steps to select **Hit** and **CompoundName** for the **Field Caption** of the cell in the first row, and **Hit** and **CompoundName** for the **Field Value** of the cell below it. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Left**.

The screenshot shows the report template editor interface. At the top, there is a header area with the placeholder text "{L:Report Title}" and the Agilent logo. Below this is a table with the following rows:

{C:DataPathName}	{V:DataPathName}
{C:DataFileName}	{V:DataFileName}
{C:SampleName}	{V:SampleName}
{C:AcqMethodFileName}	{V:AcqMethodFileName}
{C:AcqDateTime}	{V:AcqDateTime}

Below the table is a large grey area labeled "SampleChromatogram". At the bottom, there is a table with the following rows:

Peak#	RT	{C:CompoundName}							
Expression..	RetentionTir	{V:CompoundName}							

At the bottom right, there is a footer area with the placeholder text "{L:Page &[Page] of &[Pages]}" and "{L:Generated at &[Time] on &[Date]}".

- 31 Select the cell in the first row of the fourth column, and click **Textbox**.

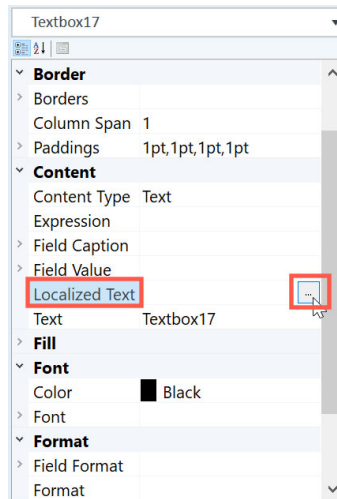
The screenshot shows the report template editor interface with the ribbon visible at the top. The ribbon has several tabs: File, Home, Font, Alignment, Colors, Page, and Image. The "Page" tab is selected, and the "Textbox" button is highlighted with a red box. Below the ribbon is the same report template editor interface as in the previous screenshot, but with a table at the bottom that has a cell selected in the first row of the fourth column.

Peak#	RT	{C:CompoundName}							
Expression..	RetentionTir	{V:CompoundName}							

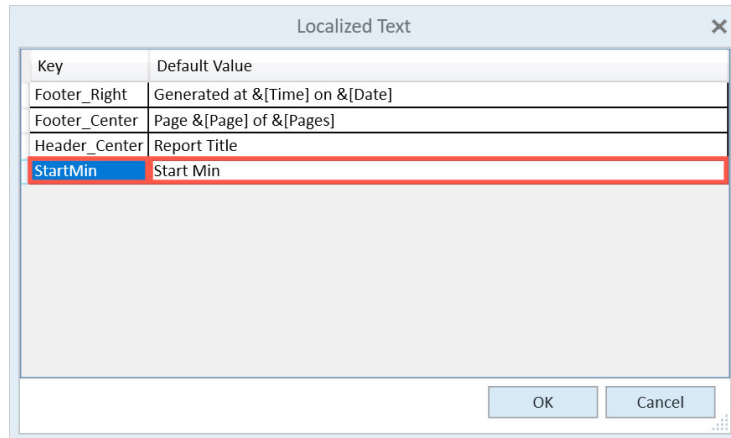
2 Report Template Workflows

Add a Compound Results table

32 Click **Localized Text** in the Properties pane, then click the ellipsis.




33 On the fourth row, enter **StartMin** in the **Key** column, and **Start Min** in the **Default Value** column. Click **OK**.




2 Report Template Workflows

Add a Compound Results table

34 In **Textbox Properties**, bold the text, and set the horizontal alignment to **Right**.

{L:Report Title}			
 Agilent <small>Trusted Answers</small>			
{C:DataPathName}	{V:DataPathName}		
{C:DataFileName}	{V:DataFileName}		
{C:SampleName}	{V:SampleName}		
{C:AcqMethodFileName}	{V:AcqMethodFileName}		
{C:AcqDateTime}	{V:AcqDateTime}		
SampleChromatogram			
Peak#	RT	{C:CompoundName}	Start Min
Expression..	RetentionTi	{V:CompoundName}	{V:StartX}
{L:Page &[Page] of &[Pages]}			{L:Generated at &[Time] on &[Date]}

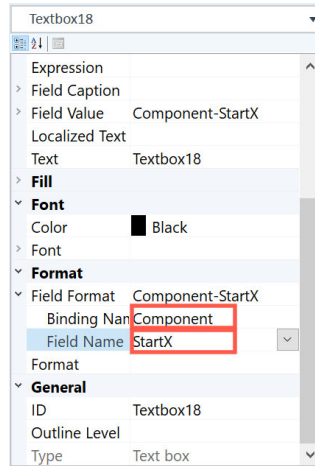
35 For the cell in the second row of the fourth column, select **Component** and **StartX** for **Field Value**, set the horizontal alignment to **Right**, and click **OK**.

{L:Report Title}			
 Agilent <small>Trusted Answers</small>			
{C:DataPathName}	{V:DataPathName}		
{C:DataFileName}	{V:DataFileName}		
{C:SampleName}	{V:SampleName}		
{C:AcqMethodFileName}	{V:AcqMethodFileName}		
{C:AcqDateTime}	{V:AcqDateTime}		
SampleChromatogram			
Peak#	RT	{C:CompoundName}	Start Min
Expression..	RetentionTi	{V:CompoundName}	{V:StartX}
{L:Page &[Page] of &[Pages]}			{L:Generated at &[Time] on &[Date]}

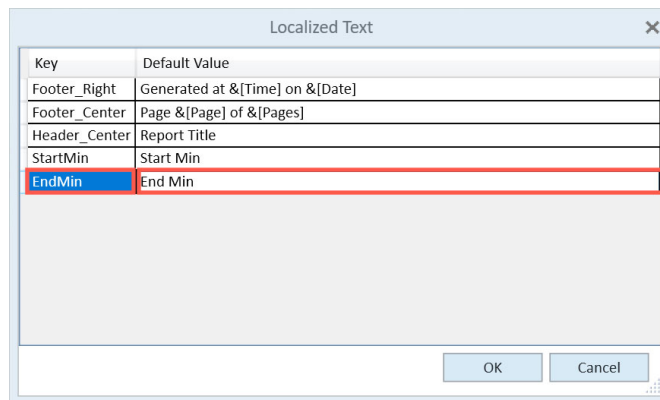
2 Report Template Workflows

Add a Compound Results table

- 36 Select the **{V:StartX}** cell, double-click **Field Format** in the Properties pane, and select **Component** from the **Binding Name** list and **StartX** from the **Field Name** list.



- 37 For the cell in the first row of the fifth column, in the **Localized Text** window, enter **EndMin** in the **Key** column, and **End Min** in the **Default Value** column, and click **OK**.



2 Report Template Workflows

Add a Compound Results table

- 38 For the cell below it, select **Component** and **EndX** for **Field Value**. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Right**.

The screenshot shows a report template with the following structure:

Peak#	RT	{C:CompoundName}	Start Min	{L:End Mii
Expression..	RetentionTiri	{V:CompoundName}	{V:StartX}	{V:EndX}

Page: {L:Page &{Page} of &{Pages}} | Generated at {L:Time} on {L:Date}

- 39 Select the cell containing **{V:EndX}**, double-click **Field Format** in the Properties pane, and select **Component** from the **Binding Name** menu, and **EndX** from the **Field Name** menu.

The screenshot shows the 'Field Format' properties pane for a field named 'Textbox20'. The 'Field Format' section is expanded, showing the following settings:

- Field Format: Component-EndX
- Binding Name: Component
- Field Name: EndX

2 Report Template Workflows

Add a Compound Results table

- 40 For the cells in the sixth column, enter **PeakArea** for **Key**, and **Peak Area** for **Default Value** in the **Localized Text** window for the cell in the first row, and select **Component** and **Area** for the **Field Value** of the cell in the second row. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Right**.

Peak#	RT	{C:CompoundName}	Start Min	{L:End Min}	Peak Area	{V:Area}			
{C:DataPathName}	{V:DataPathName}	{C:DataFileName}	{V:DataFileName}	{C:SampleName}	{V:SampleName}	{C:AcqMethodFileName}	{V:AcqMethodFileName}	{C:AcqDateTime}	{V:AcqDateTime}

- 41 Select the **{V:Area}** cell, and select **Component** and **Area** in the **Field format** of the Properties pane.

Textbox22

> Paddings 1pt,1pt,1pt,1pt

> Content

Content Type FieldValue

Expression

> Field Caption

> Field Value Component-Area

Localized Text

Text Textbox22

> Fill

> Font

Color Black

> Font

> Format

> Field Format Component-Area

Binding Name Component

Field Name Area

Format

> General

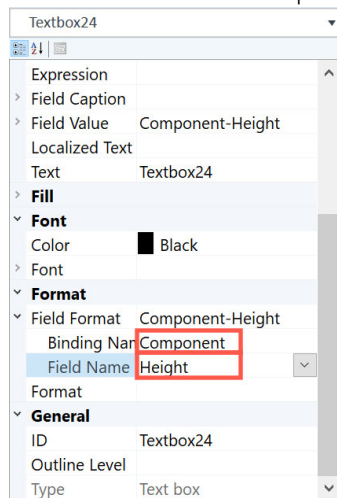
2 Report Template Workflows

Add a Compound Results table

- 42 For the cells in the seventh column, enter **PeakHeight** for **Key**, and **Peak Height** for **Default Value** in the **Localized Text** window for the cell in the first row, and select **Component** and **Height** for the **Field Value** of the cell in the second row. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Right**.

{L:Report Title}						
{C:DataPathName}	{V:DataPathName}					
{C:DataFileName}	{V:DataFileName}					
{C:SampleName}	{V:SampleName}					
{C:AcqMethodFileName}	{V:AcqMethodFileName}					
{C:AcqDateTime}	{V:AcqDateTime}					
SampleChromatogram						
Peak#	RT	{C:CompoundName}	Start Min	{L:End Min}	Peak Area	Peak Height
Expression...	RetentionTime	{V:CompoundName}	{V:StartX}	{V:EndX}	{V:Area}	{V:Height}
			{L:Page &[Page] of &[Pages]}		{L:Generated at &[Time] on &[Date]}	

- 43 Select the **{V:Height}** cell, and select **Component** and **Height** in the **Field format** section of the Properties pane.



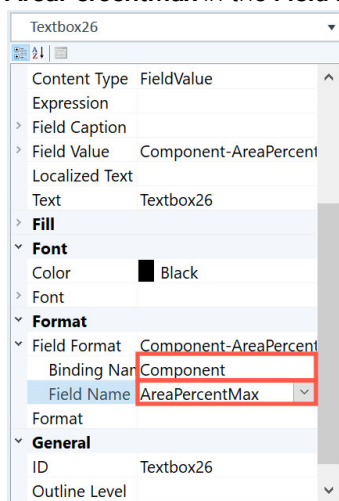
2 Report Template Workflows

Add a Compound Results table

- 44 For the cells in the eighth column, enter **Peak%Max** for **Key**, and **Peak % Max** for **Default Value** in the **Localized Text** window for the cell in the first row, and select **Component** and **AreaPercentMax** for the **Field Value** of the cell in the second row. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Right**.

		{L:Report Title}		Agilent	Trusted Analysis
{C:DataPathName}		{V:DataPathName}			
{C:DataFileName}		{V:DataFileName}			
{C:SampleName}		{V:SampleName}			
{C:AcqMethodFileName}		{V:AcqMethodFileName}			
{C:AcqDateTime}		{V:AcqDateTime}			
SampleChromatogram					
Peak#	RT	{C:CompoundName}	Start Min	{L:End Min	Peak Area
Expression...	RetentionTime	{V:CompoundName}	{V:StartX}	{V:EndX}	{V:Area}
					{V:Height} ercentMax}
				{L:Page &[Page] of &[Pages]}	{L:Generated at &[Time] on &[Date]}

- 45 Select the **{V:AreaPercentMax}** cell, and select **Component** and **AreaPercentMax** in the **Field format** section of the Properties pane.



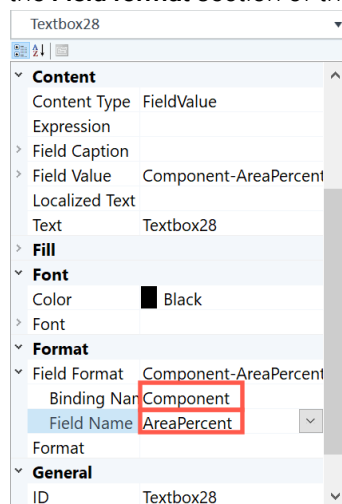
2 Report Template Workflows

Add a Compound Results table

- 46 For the cells in the ninth column, enter **Peak%Total** for **Key**, and **Peak % Total** for **Default Value** in the **Localized Text** window for the cell in the first row, and select **Component** and **AreaPercent** for the **Field Value** of the cell in the second row. Bold the text of the cell in the first row, and change the horizontal alignment of both cells to **Right**.

		{L:Report Title}		Agilent <small>Trusted Accounts</small>
{C:DataPathName}		{V:DataPathName}		
{C:DataFileName}		{V:DataFileName}		
{C:SampleName}		{V:SampleName}		
{C:AcqMethodFileName}		{V:AcqMethodFileName}		
{C:AcqDateTime}		{V:AcqDateTime}		
SampleChromatogram				
Peak#	RT	{C:CompoundName}	Start Min	{L:End Min Peak Area} k % Max
Expression..	RetentionTime	{V:CompoundName}	{V:StartX} {V:EndX}	{V:Area} {V:Height} ercentMax} reaPercent}
			{L:Page &{Page} of &{Pages}}	{L:Generated at &{Time} on &{Date}}

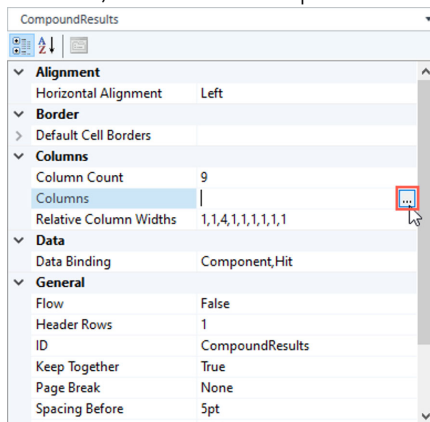
- 47 Select the **{V:AreaPercent}** cell, and select **Component** and **AreaPercent** in the **Field format** section of the Properties pane.



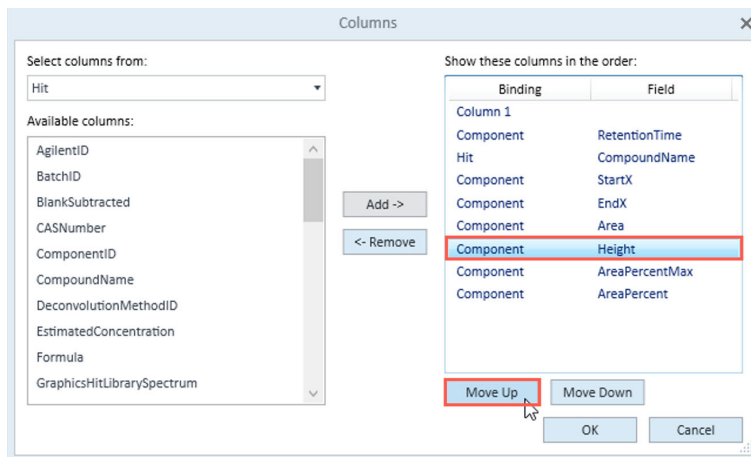
2 Report Template Workflows

Add a Compound Results table

- 48 Select **CompoundResults** from the list at the top of the Properties pane. Click **Columns**, then click the ellipsis.



- 49 In the **Columns** window, select **Height**, and click **Move Up**. Click **OK**.

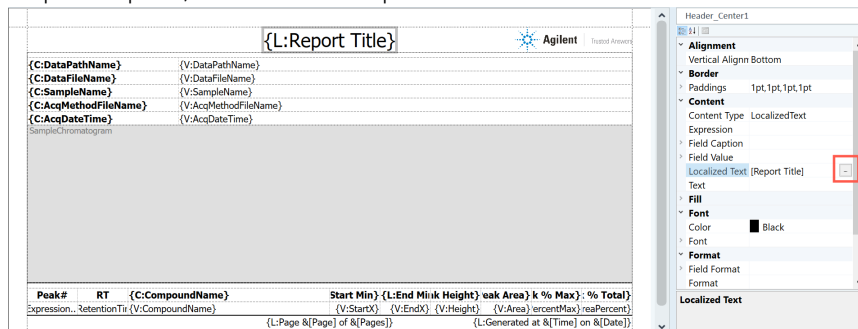


2 Report Template Workflows

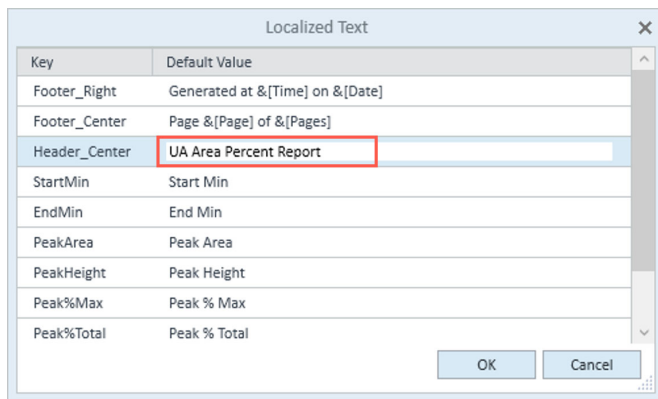
Add a report title

Add a report title

- 1 Select the **{L:Report Title}** cell in the page header. Click **Localized Text** in the Properties pane, then click the ellipsis.



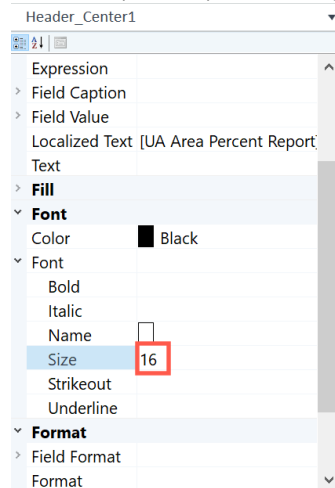
- 2 Enter **UA Area Percent Report** in the **Default Value** column for **Report Title**. Click **OK**.



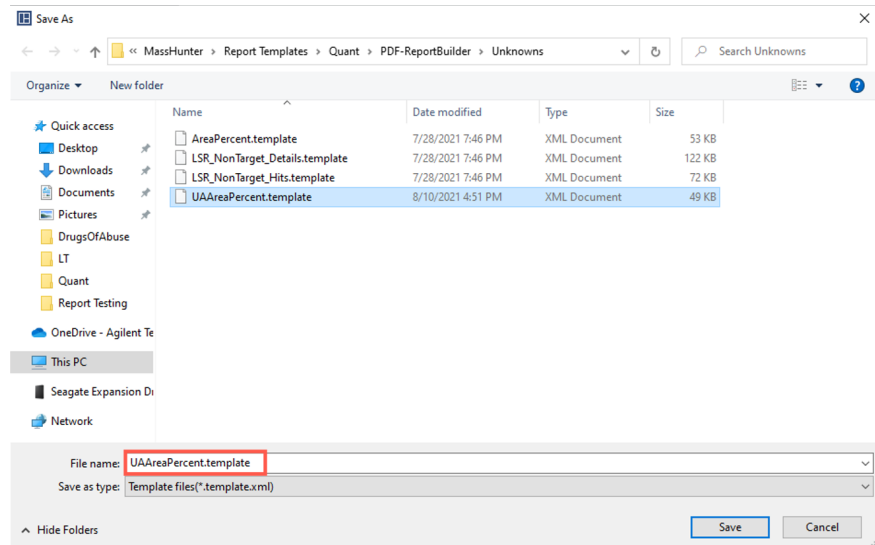
2 Report Template Workflows

Add a report title

- 3 In the Properties pane, change the font size to **16**.



- 4 Click **File > Save As**, and save the template as **UAAreaPercent.template.xml**.



2 Report Template Workflows

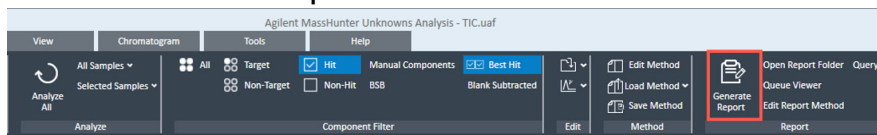
Generate a report

Generate a report

- 1 Open a TIC analysis in MassHunter Unknowns Analysis.



- 2 Click **Home > Generate Report**.



2 Report Template Workflows

Generate a report

- 3 Under **Report method**, click **Choose**, navigate to the relevant folder, and load your method.

Report

Report folder: _____
X:\Sue Report Testing\TestingData\RI-PEST-MATRIX\Unknowns Browse...

Report method: _____
X:\Sue Report Testing\TestingData\RI-PEST-MATRIX\AreaPercent.m
Choose... New... Edit...

Samples: _____
 All samples
 Selected sample(s)

Generate: _____
 Generate reports now
 Open report folder after reports generated
 Queue report task
 Start Queue Viewer

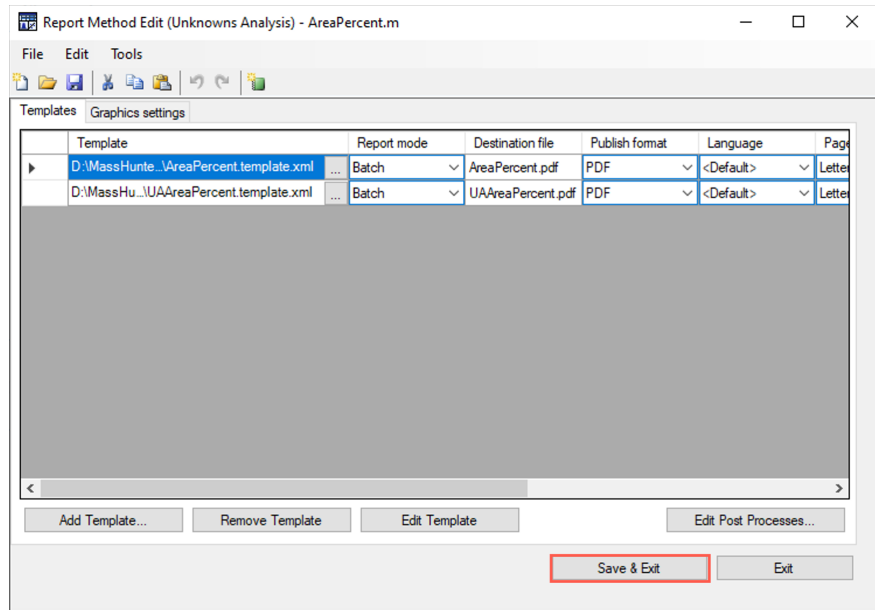
OK Cancel

- 4 Under **Report method**, click **Edit**.
- 5 Click **Add Template**.
- 6 Search for and select your report template, and click **Open**.

2 Report Template Workflows

Generate a report

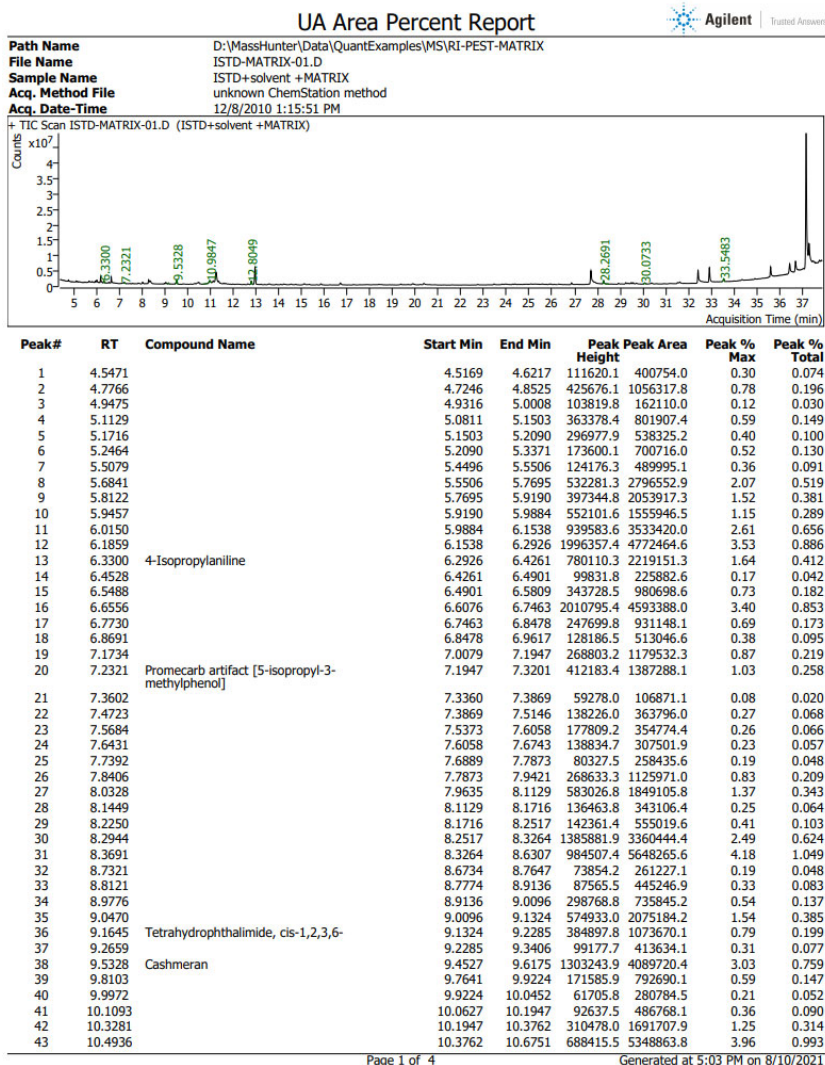
7 Click **Save & Exit**.



2 Report Template Workflows

Generate a report

- Click **OK** at the bottom of the **Report** window. A PDF report is generated using the selected batch files.



3 Modify a Report Template

- Replace the Agilent Logo with Your Own 98
- Replace a Column in the Report 99
- Replace a column using the Textbox Properties window 101
- Remove a Column From the Report 103
- Adjust the Width of a Column 104
- Filter Results Using Special Conditions 105
- Adjust the Size of a Graphic 107
- Add a New Column to an Existing Report 108
- Edit an Expression 112
- Preview the Report 113

In this section you will modify a report template by completing a series of tasks, such as replacing the logo, modifying tables, filtering results, adjusting graphic sizes, and adding an expression.

3 Modify a Report Template

Replace the Agilent Logo with Your Own

Replace the Agilent Logo with Your Own

- 1 Open the **Report Builder**.
- 2 Select **File > Open**.
- 3 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 4 The Agilent logo is in the upper right corner of the report template. This is the header section of the report page.

Select the logo graphic. The filename **AgilentLogo.tif** appears in the right panel.
- 5 If the logo you want to use is in the same file location as the report template, enter the file name in the **File** field, and press **Enter**.

If the logo you want to use is in a different location than the report template, enter the full path of the file in the **File** field, and press **Enter**.
- 6 Save the template.

3 Modify a Report Template

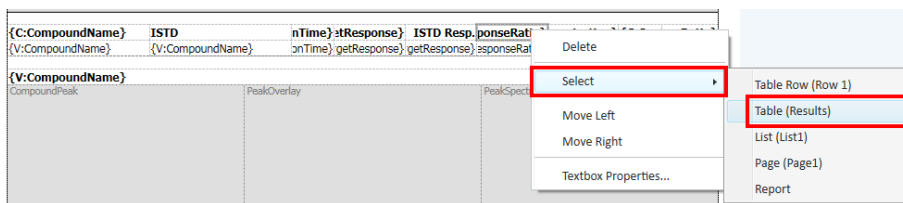
Replace a Column in the Report

Replace a Column in the Report

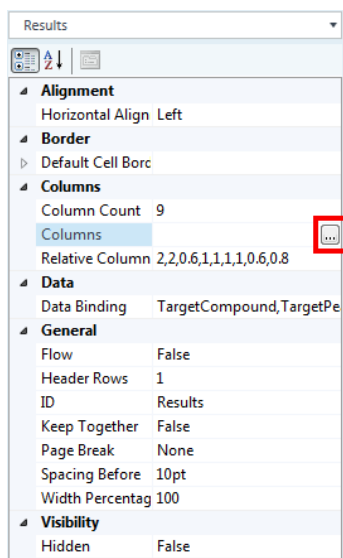
You can replace a column in a table through either the **Columns** window or the **Textbox Properties** window.

Replace a column using the Columns window

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Right-click the **ResponseRatio** column header, and select **Select > Table (Results)**.



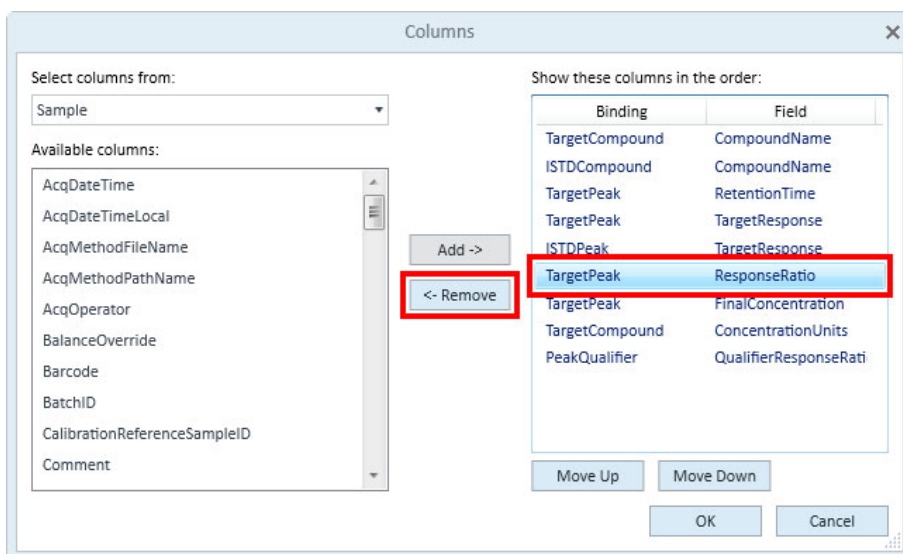
- 3 In the **Columns** field in the **Properties** pane, click the ellipsis.



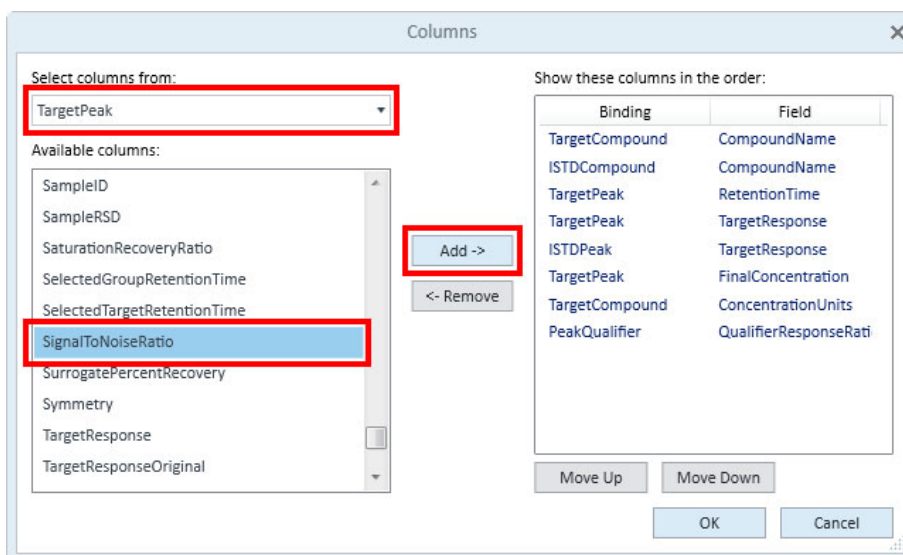
3 Modify a Report Template

Replace a column using the Columns window

- 4 In the **Columns** window, select **ResponseRatio**, and click **Remove**.



- 5 Select **TargetPeak** from the **Select columns from:** list, select **SignalToNoiseRatio** in the **Available columns** pane, and click **Add**.



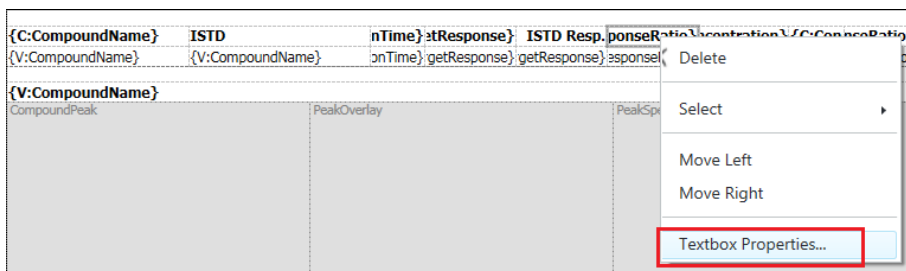
- 6 Click **Move Up** or **Move Down** to move the column to the desired location.
- 7 Click **OK**. The **ResponseRatio** column is replaced by the **SignalToNoiseRatio** column.

3 Modify a Report Template

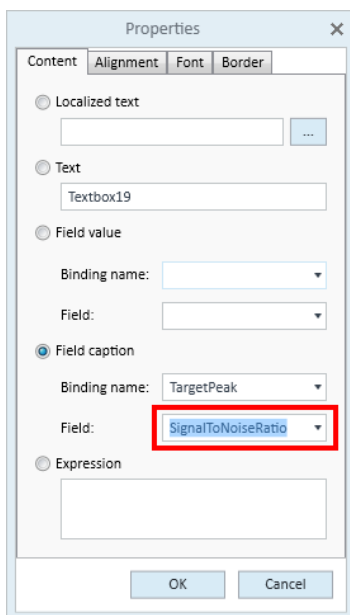
Replace a column using the Textbox Properties window

Replace a column using the Textbox Properties window

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Right-click the **ResponseRatio** column header, and select **Textbox Properties**.



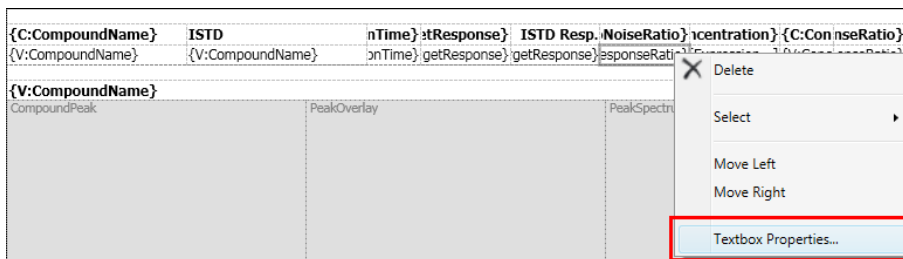
- 3 Select **SignalToNoiseRatio** from the **Field Caption > Field** list, and click **OK**.



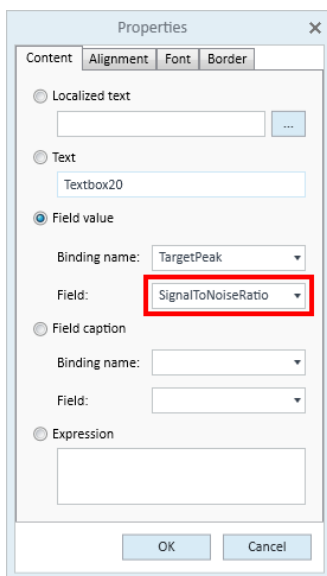
3 Modify a Report Template

Replace a column using the Textbox Properties window

4 Right-click the **ResponseRatio** value, and select **Textbox Properties**.



5 Select **SignalToNoiseRatio** from the **Field Value > Field** list, and click **OK**



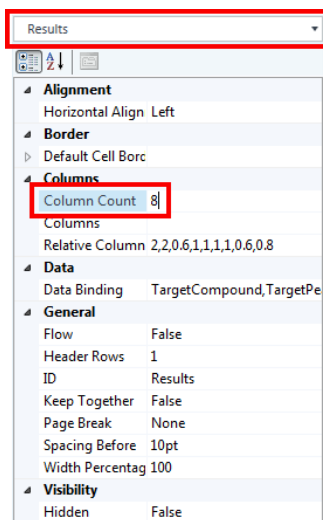
6 The **ResponseRatio** column is replaced by the **SignalToNoiseRatio** column. Save the template.

3 Modify a Report Template

Remove a Column From the Report

Remove a Column From the Report

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Select the table that contains the **QualifierResponseRatio** column (Results).
- 3 In the **Column Count** field in the **Properties** pane, change the column count from 9 to **8**.



The last column in the table, **QualifierResponseRatio**, is removed.

- 4 Save the template.

To remove a column that is not the last column in the table:

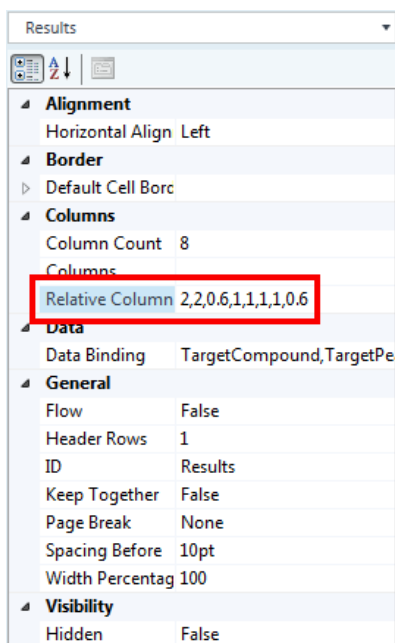
- 1 Right-click the column to remove, and select **Move Right**.
Repeat this step until the column is the last column in the table.
- 2 In the **Column Count** field in the **Properties** pane, reduce the column count by 1 to remove the last column.
- 3 Save the template.

3 Modify a Report Template

Adjust the Width of a Column

Adjust the Width of a Column

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Select a table.
- 3 In the **Relative Column Widths** field in the **Properties** pane, change the values to adjust the widths of the columns in the table.



- 4 Save the template.

3 Modify a Report Template

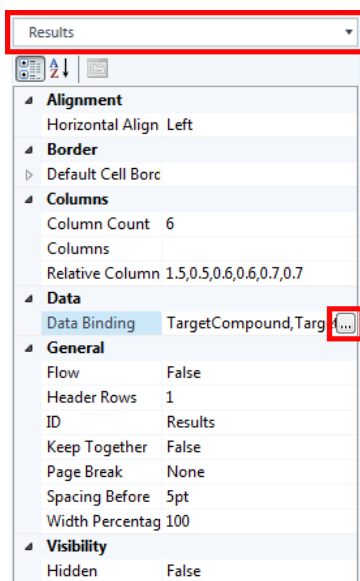
Filter Results Using Special Conditions

Filter Results Using Special Conditions

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Sample_AboveLOQ.template.xml**.

This report filters results lower than the LOQ. Use this procedure to modify the data binding to filter results lower than the LOD.

- 2 Select the **Results** table in the **Properties** pane, and click the ellipsis in the **Data Binding** field.
- 3 In the **Data Binding** window, select **Target Peak**, and then select **AND <CalculatedConcentration> >= <TargetCompound:LimitOfQuantitation>**.



3 Modify a Report Template

Filter Results Using Special Conditions

4 Click **Edit**.

The screenshot shows the 'Data Binding' dialog box. On the left, a list of data bindings includes 'TargetCompound' and 'TargetPeak', with 'TargetPeak' selected. The 'Binding name' is 'TargetPeak' and the 'Data name' is 'Peak'. The 'Filters' section contains several conditions, with the last one, 'AND <CalculatedConcentration> >= <TargetCompound.LimitOfQuantitation>', highlighted. The 'Orders' section is empty. The 'Expression' field is also empty. Buttons for 'Add', 'Delete', 'Move Up', 'Move Down', 'New...', 'Edit...', and 'Delete' are visible. The 'Left outer join' checkbox is unchecked. 'OK' and 'Cancel' buttons are at the bottom right.

5 Select **LimitOfDetection** from the **Field name** list, and click **OK**.

The screenshot shows the 'Data Binding Filter' dialog box. The 'Field' is 'CalculatedConcentration', the 'Operator' is '>=', and the 'Concatenation type' is 'AND'. The 'Fixed value' field is empty. The 'Bound value' field is empty. The 'Binding name' is 'TargetCompound'. The 'Field name' is 'LimitOfDetection', which is highlighted. 'OK' and 'Cancel' buttons are at the bottom.

6 Click **OK**.

7 Save the template.

3 Modify a Report Template

Adjust the Size of a Graphic

Adjust the Size of a Graphic

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Select the graphic to resize.
- 3 In the **Height** and **Width Percentage** fields in the **Properties** pane, change the values to adjust the size of the graphic.

The screenshot displays the Report Builder interface. On the left, the XML template is shown with the following content:

```
{C:AcqDateTime} {V:AcqDateTime} {C:DataFileName} {V:Dat
{C:SampleType} {V:SampleType} {C:SampleName} {V:Sam
{C:Dilution} {V:Dilution} {C:AcqMethodFileName} {V:Acc

{L:Sample Chromatogram}
SampleChromatogram

{C:CompoundName} ISTD nTime}stResponse} ISTD Resp.ponseRatio} cent
{V:CompoundName} {V:CompoundName} onTime}.getResponse}.getResponse}.sponseRatio}. Expre
{V:CompoundName}
CompoundPeak PeakOverlay PeakSpectrum
```

On the right, the Properties pane is open, showing the following settings:

- General**
 - Height: 180pt
 - ID: SampleChromatogram1
 - Name: SampleChromatogram
 - Width Percentage: 100
- ID**
 - BatchID: Parameter [BatchID]
 - SampleID: Parameter [SampleID]
- Misc**
 - Borders: [empty]
 - Column Span: 1
- MS**
 - MSScanType: [empty]
- Signal**
 - DeviceName: Parameter [DeviceName]
 - OrdinalNumber: Parameter [OrdinalNumber]
 - SignalName: Parameter [SignalName]

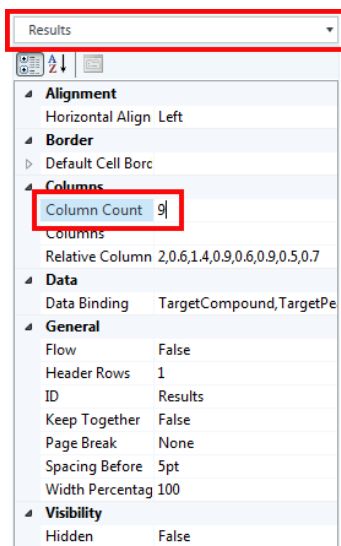
- 4 Save the template.

3 Modify a Report Template

Add a New Column to an Existing Report

Add a New Column to an Existing Report

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ESTD.template.xml**.
- 2 Select the **Results** table from the **Properties** pane, and change the **Column Count** value from 8 to **9**.



A new empty column is added to the end of the table.

3 Modify a Report Template

Add a New Column to an Existing Report

- 3 Add a textbox to the empty column header and value. Separately select the empty cells in the table, and click **Textbox**.

The screenshot shows the 'Home' ribbon of the Report Builder software. The 'Textbox' button is highlighted with a red box. Below the ribbon, a table contains various report fields. A red line points from the 'Textbox' button to the table area.

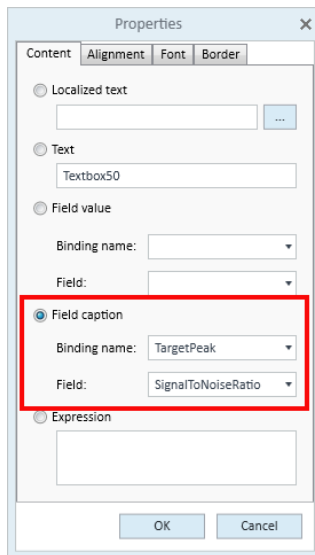
{C:AnalysisTimeStamp}	{V:AnalysisTimeStamp}	{C:AnalystName}	{V:AnalystName}
{L:Report Generation Time}	&[DateTime]	{C:ReportGeneratorName}	[Expression...]
{L:Calibration Last Update}	{V:CalibrationLastUpdatedTimeSta	{C:BatchState}	{V:BatchState}
{C:AnalyzeQuantVersion}	{V:AnalyzeQuantVersion}	{C:ReportQuantVersion}	&[CurrentVersion]
{C:AcqDateTime}	{V:AcqDateTime}	{C:DataFileName}	{V:DataFileName}
{C:SampleType}	{V:SampleType}	{C:SampleName}	{V:SampleName}
{C:Dilution}	{V:Dilution}	{C:AcqMethodFileName}	{V:AcqMethodFileName}
{L:Sample Chromatogram}			
SampleChromatogram			
{C:CompoundName}	{C:Reten	{C:Transition}	{C:TargetRe: {C:Height}ncentration} {C:ConcnseRatio} Textbox50
{V:CompoundName}	{V:Retent	{V:Transition}	{V:TargetResp {V:Height [Expression...]} {V:ConcnseRatio}

- 4 Right-click the column header textbox and select **Textbox Properties**.

3 Modify a Report Template

Add a New Column to an Existing Report

- 5 Select **Field caption**, select **TargetPeak** from the **Binding name** list, and then select **SignalToNoiseRatio** from the **Field** list.



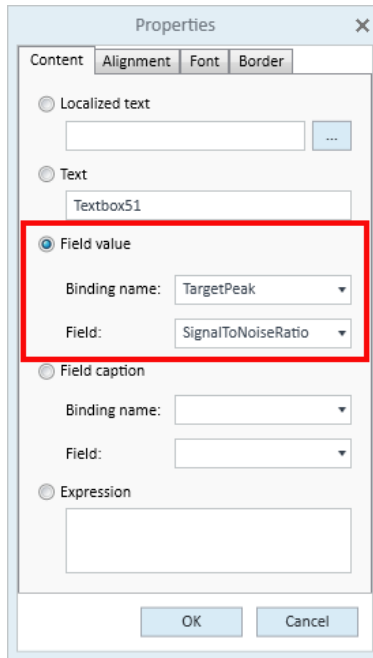
The image shows a 'Properties' dialog box with a close button (X) in the top right corner. It has four tabs: 'Content', 'Alignment', 'Font', and 'Border'. The 'Content' tab is active. There are five radio button options: 'Localized text', 'Text', 'Field value', 'Field caption', and 'Expression'. The 'Field caption' option is selected and highlighted with a red rectangular box. Below this option, there are two dropdown menus: 'Binding name:' which is set to 'TargetPeak', and 'Field:' which is set to 'SignalToNoiseRatio'. At the bottom of the dialog are 'OK' and 'Cancel' buttons.

- 6 Click **OK**.
- 7 Right-click the column value textbox and select **Textbox Properties**.

3 Modify a Report Template

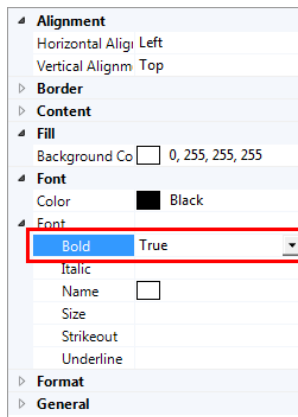
Add a New Column to an Existing Report

- 8 Select **Field value**, select **TargetPeak** from the **Binding name** list, and then select **SignalToNoiseRatio** from the **Field** list.



The screenshot shows the 'Properties' dialog box with the 'Field value' section selected. The 'Binding name' dropdown is set to 'TargetPeak' and the 'Field' dropdown is set to 'SignalToNoiseRatio'. The 'Field value' section is highlighted with a red box.

- 9 Select **True** for the **Bold** field in the **Properties** pane for the field caption and field value.



The screenshot shows the 'Properties' dialog box with the 'Font' section selected. The 'Bold' field is set to 'True' and is highlighted with a red box.

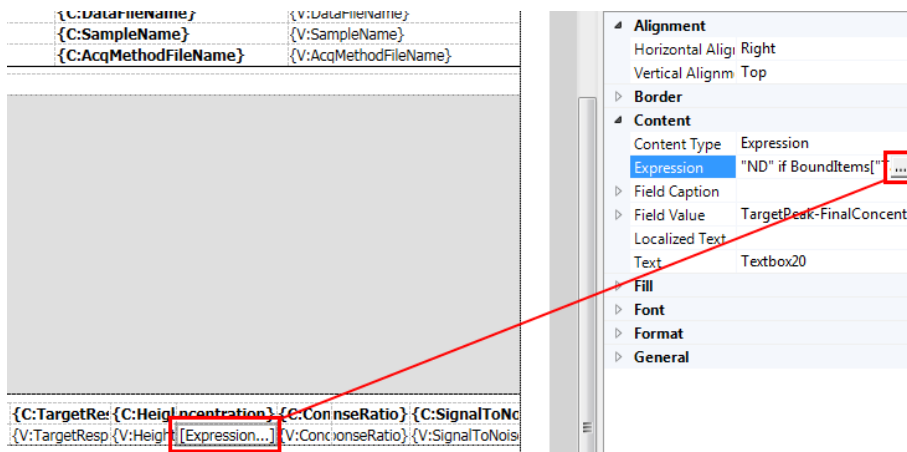
- 10 Save the template.

3 Modify a Report Template

Edit an Expression

Edit an Expression

- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Select the expression in the **FinalConc** column in **Table 2**, and click the ellipsis for the **Expression** field in the **Properties** pane.



- 3 Modify the expression as needed in the **ExpressionWindow** window, and click **OK**.
- 4 Save the template.

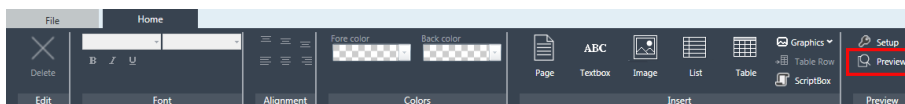
3 Modify a Report Template

Preview the Report

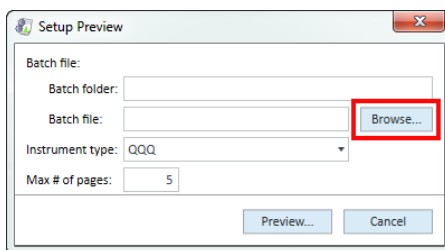
Preview the Report

After changes have been made to the report template, you can preview the report by setting up the Report Builder with your dataset.

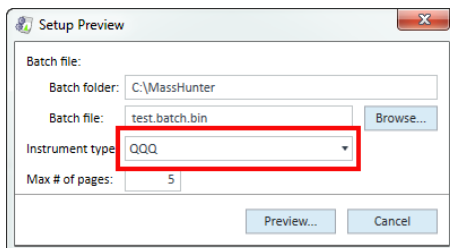
- 1 Navigate to **D:\MassHunter\Report Templates\Quant\PDF-ReportBuilder\Quant**, and open **Gen_Samples_ISTD.template.xml**.
- 2 Click **Preview**.



- 3 In the **Setup Preview** window, click **Browse**, then navigate to and select your batch file.



- 4 Select the **Instrument type** to match your dataset.

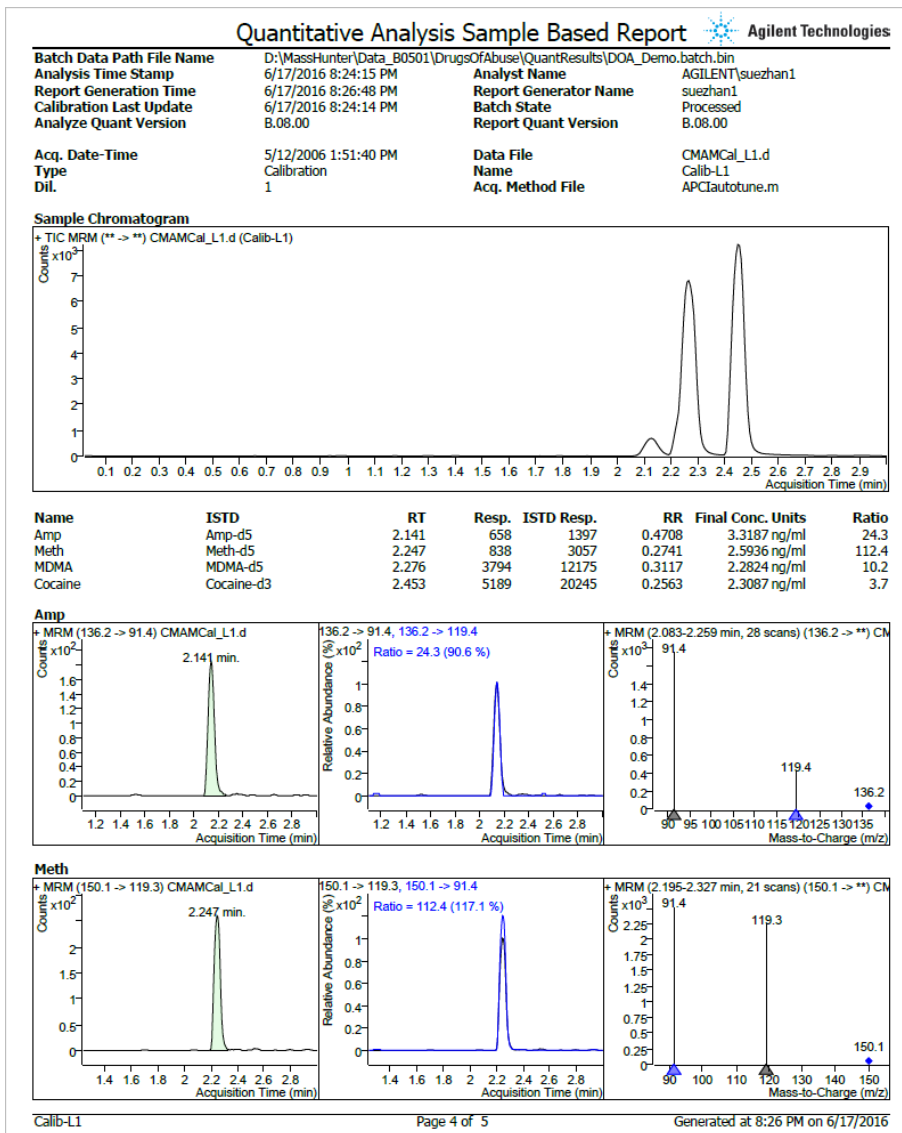


- 5 Click **Preview**. A PDF report is generated using the batch file selected.

3 Modify a Report Template

Preview the Report

Navigate the report page by page. Below is an example of page 4 using the **DrugsofAbuse** dataset.



www.agilent.com

© Agilent Technologies, Inc. 2021

November 2021



G3336-90061

