



Agilent MassHunter Drug Analysis Mode Using Quantitative Analysis

Workflow Guide



Review the Intelligent Sequencing Feature

Create the Data Acquisition Method

Create the Quant Analysis Method

Run Samples



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2013

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

G6845-90007

Edition

First edition, August 2013

Printed in USA

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

Warranty

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

Technology Licenses

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

Restricted Rights Legend

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

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

Safety Notices

CAUTION

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

WARNING

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

In This Guide...

This workflow guide describes how a method can be created to perform data acquisition in MassHunter GCMS Acquisition and data analysis in MassHunter Quantitative Analysis, using the Drug Analysis Workflow mode. Also, for those users interested in migrating from MSD ChemStation to MassHunter Quantitative Analysis, Chapter 4 includes a section on converting an existing MSD ChemStation Method to a MassHunter Quantitative Analysis method.

This document emphasizes the use and understanding of Intelligent Sequencing features in MassHunter Drug Workflow mode, which are not currently covered in other documents created for MassHunter GCMS Acquisition and Quantitative Analysis.

More common operations, not directly associated with Drug Analysis Workflow mode and Intelligent Sequencing, such as setting GC Parameters and real-time plot displays, are briefly discussed here, but are covered in more detail in both online Help and Familiarization Guides. Please refer to the online Help for more details on these topics and for a link to an unabridged version of the MassHunter GCMS Acquisition Software for 5975/5977 Series GC/MSD Familiarization Guide (G1701-90110).

A brief summary of chapter contents for this Workflow Guide follows.

1 Before You Begin

Chapter 1 describes how to set up your MassHunter GCMS Acquisition and MassHunter Quantitative Data Analysis programs for using the Drug Workflow mode user interface (UI). Items specific to this UI are then reviewed.

2 About Intelligent Sequencing

Chapter 2 describes Intelligent Sequencing and how it uses keywords to initiate actions based on results from Data Analysis. These actions include injecting blanks to reduce sample carryover, reinjecting a specimen including reducing the injection volume when its concentration is over a predefined limit, skipping to another batch, or just waiting for an operator to intervene. Because a number of actions may be employed within a given sequence, flow diagrams are included to aid in following the decision process.

3 Create the Data Acquisition Method

Chapter 3 describes how to set up a SIM method for data acquisition. A Data Acquisition method must exist prior to the creation of a Quantitative Data Analysis method that uses Intelligent Sequencing.

4 Create the Quantitative Analysis Method

Chapter 4 describes how to create a Drug Quant method, compatible with Intelligent Sequencing, by editing an existing method. It starts out by describing how you can translate an existing MSD ChemStation method into a MassHunter Quantitative Analysis method, if desired. It then proceeds with setting up calibration curves for

5 Run Samples

various drugs and setting parameters required by Intelligent Sequencing. This includes setting the allowable specifications used during analysis and the set up of control samples used for monitoring the calibration state.

6 Update the Calibration

Chapter 5 explains how to set up and run a Sequence Table with multiple batches of samples using the keyword **NewBatch**. This keyword indicates where MassHunter should continue processing when a *Skip to next batch* command is generated by Intelligent Sequencing.

Chapter 6 describes how the Sequence Table Editor (STE) in MassHunter GCMS Acquisition can be used to automate the process of updating the calibration table. As required by governing regulations, the calibration stored in a method must be updated when a specified time has elapsed or when a quality control sample indicates an unacceptable deviation from the stored calibration curve.

Where to Find More Information

Accompanying your hardware and software is a comprehensive collection of manuals, videos, user applications, and method development tools. These are located on the:

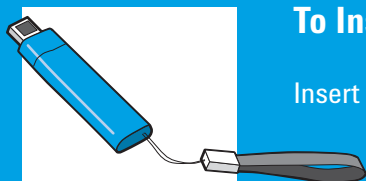
- Agilent GC and GC/MS Manuals and Tools DVD set
- Agilent GC/MS Software Information and Manuals memory stick



To Install Your Hardware Library

Insert Disk 1 into your DVD drive and follow the prompts.

This can be installed by anyone who has authority to copy information onto the receiving computer.



To Install Your Software Library

Insert the memory stick into a USB port and follow the prompts.

This can be installed by anyone who has authority to copy information onto the receiving computer.

Contents

1 Before You Begin

- Configure MassHunter GCMS Acquisition for Drug Analysis 8
- Configure MassHunter Quant for Drug Analysis 10
- Understand the Directory Structure 12
- Review the Batch Menu 13

2 About Intelligent Sequencing

- What is Intelligent Sequencing? 16
- How does it work? 16
- Where are limits and parameters defined? 16
- What actions can be taken by Intelligent Sequencing? 18
- What are some examples of how Intelligent Sequencing can be used? 19
- How are Blanks handled? 20
- How are Negatives and Controls handled? 24
- How are Specimens handled? 30
- Maximum # of Retries 37

3 Create the Data Acquisition Method

- Step 1: Load a data acquisition method 40
- Step 2: Select the parts of the method to edit 40
- Step 3: Describe the method and where it is saved 41
- Step 4: Complete the Intelligent Sequencing parameters 42
- Step 5: Complete the Standard Instrument Acquisition dialogs 43
- Step 6: Create a SIM method from a Scan method 44
- Step 7: Save the Method 45

4 Create the Quantitative Analysis Method

- Introduction 48
- Step 1: Convert an MSD ChemStation method (Optional) 48
- Step 2: Set up the method for Drug Quant 50
- Step 3: Change each compound's Type from Scan to SIM 53
- Step 4: Add a multi-level calibration to the method 54
- Step 5: Add Control Samples to the Method 57
- Step 6: Add ISTD Compounds to Sample as Target Compounds 59
- Step 7: Set up outliers for intelligent sequencing decisions 61
- Step 8: Save and Validate the Method 67

5 Run Samples

Introduction 70

Step 1: Load the default Sequence 70

Step 2: Edit the Sequence Table 71

Step 3: Specify reports 73

Step 4: Save the Sequence Table 73

Step 5: Run the Sequence 74

6 Update the Calibration

Introduction 76

Step 1: Create the Sequence Table 76

Step 2: Prepare the calibration samples 80

Step 3: Prepare the cutoff calibration sample 80

Step 4: Load the calibration sample vials in the ALS 80

Step 5: Run the Calibration Sequence 81

1 Before You Begin

Configure MassHunter GCMS Acquisition for Drug Analysis 8

Configure MassHunter Quant for Drug Analysis 10

Understand the Directory Structure 12

Review the Batch Menu 13

Before You Begin

Configure Drug
Analysis

Review Items
Specific to the
Drug Analysis
Mode



Agilent Technologies

Configure MassHunter GCMS Acquisition for Drug Analysis

1. Double-click the GCMS Configuration desktop icon to launch the Agilent GC/MS Configuration program.
2. Select the instrument name that you will be running to acquire the data. Instrument 1 is selected in this example.
3. Select the **Drug Analysis Workflow Mode** and click **OK** to close the dialog.

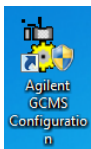
Click **Yes** to confirm the configuration and exit the Agilent GC/MS Configuration program.

Depending on your instrument, MassHunter GCMS Acquisition and MassHunter Quantitative Analysis may be set up to run in several Workflow Modes, including:

- Enhanced
- Drug Analysis
- EnviroQuant (EPA)
- Aromatics in Gasoline

Here we are going to be using the **Drug Analysis Workflow Mode**. So, before doing anything else, you must set up the MassHunter GCMS Acquisition program and the MassHunter Quantitative Analysis program to run in the Drug Analysis Workflow Mode.

To reconfigure an existing GC/MSD instrument to work in the Drug Analysis Mode:



Agilent GC/MS Instrument Configuration

File Configure Help

1 2 3 4 ?

Execute

Current Agilent GC/MS Instrument Configuration

	Name	Offline	MS	MS IP	Available Sources	GC	GC IP	Workflow Mode
1	Kermit	<input checked="" type="checkbox"/>	5977	192.168.1.201		7890	192.168.1.203	Drug Analysis
2	<none>	<input type="checkbox"/>	<none>			<none>		None
3	<none>	<input type="checkbox"/>	<none>			<none>		None
4	<none>	<input type="checkbox"/>	<none>			<none>		None

Agilent GC/MS Instrument Configuration

Instrument Name: Kermit

Laboratory ID Number: 201

Offline Instrument

Mass Spectrometer

Model: 5977

Address: 192.168.1.201

DC Polarity: Positive (+) Negative (-)

Gas Chromatograph

Model: 7890

Address: 192.168.1.203

Headspace Type: <none>

Workflow Mode: Drug Analysis

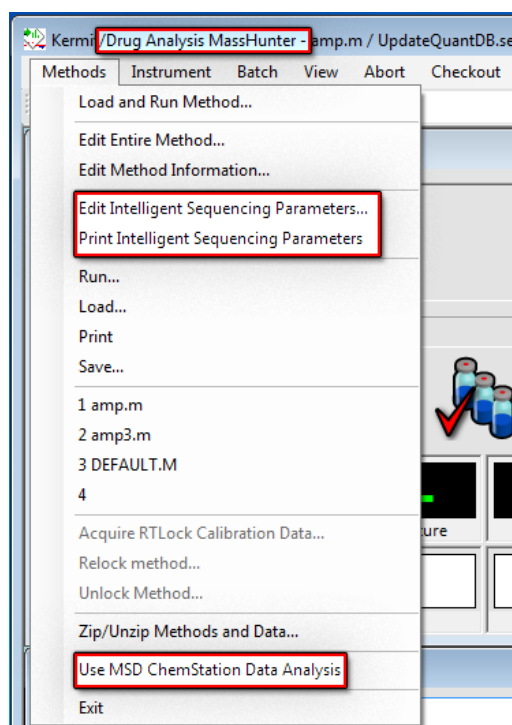
OK Cancel Help

4. Double-click the Instrument icon to launch MassHunter GCMS Acquisition.



5. Notice the items that are unique to the Drug Analysis Workflow mode.

The next time you start MassHunter GCMS Acquisition you will see Drug Analysis in the Title Bar, and under the Methods menu, you will see options for **Intelligent Sequencing Parameters**, and possibly one for **Use MSD ChemStation Data Analysis**.



Configure MassHunter Quant for Drug Analysis

Check for the Startup icon

When MassHunter Quantitative Analysis is installed, a group of icons for starting Quantitative Analysis, similar to those shown here, is placed on the windows desktop.

To begin MassHunter Quantitative Analysis you simply double-click on the applicable icon.

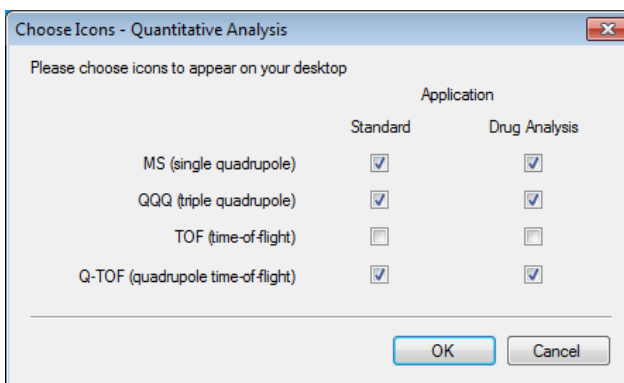


For example, to start a Quantitative Analysis session for Triple Quadrupole Data in the Drug Quant workflow mode you would click the desktop icon labeled **Drug Quant (QQQ)**. The Quant program is then optimized for Triple Quadrupole Data in the Drug Quant workflow mode.

Add a startup icon

1. From the windows Start menu select **Agilent\MassHunter Workstation\Quant Tools\Setup Desktop Icons**.
2. Check the **Drug Analysis** mode for your instrument(s).
3. Click **OK** to close the dialog and add your newly selected startup icon(s) to the windows desktop.

If you do not see a desktop icon labeled **Drug Quant** for your instrument, add it as follows.



In this example, both the Standard and Drug Analysis modes are selected for MS, Triple Quadrupole, and QTOF instruments.

Confirm the Data Analysis program can communicate with Data Acquisition program

This workflow uses:

- MassHunter GCMS Acquisition to acquire data
- MassHunter Quantitative Analysis (with Intelligent Sequencing) for Data Analysis

Intelligent sequencing uses automated processing and feedback between these two programs as it processes each sample. Therefore, be sure that:

MassHunter Quantitative Analysis is installed on the same PC as the MassHunter GCMS Acquisition program that will acquire the data.

If MassHunter Quant is installed on the same PC as the MassHunter GCMS Acquisition program, by default, it is generally configured to be used for data analysis of MassHunter GCMS Acquisition files. However, to be sure, in MassHunter GCMS Acquisition, select **Method** and, if the **Use MSD ChemStation Data Analysis** option is listed in the menu, be sure that it is **not** checked. (This menu item only exists when both MassHunter Quant and MSD ChemStation Data Analysis are installed on the same computer.)

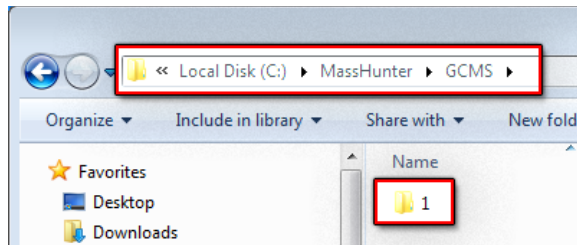
Remember, if this menu item is selected, the system will use MSD ChemStation for data analysis, and for this workflow that is not what you want. This workflow uses MassHunter Quantitative Analysis for Data Analysis.

Understand the Directory Structure

1. Locate the instrument directories.

You can configure and run up to four instruments with MassHunter GCMS Acquisition.

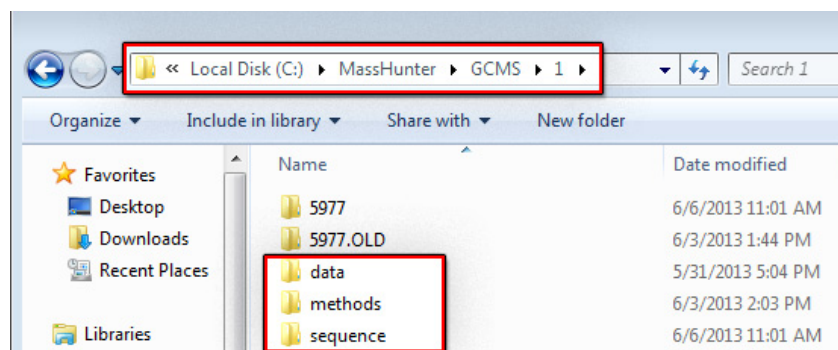
For each instrument you configure, MassHunter GCMS Acquisition will create a numbered directory corresponding to the instrument number (**drive:\MassHunter\GCMS\1** for example).



Under each instrument directory (**1** shown here), you will see a default data, methods, and sequence subdirectory, as shown in the next example.

2. Review the default data, methods, and sequence directories.

These are the recommended and default locations for your data, methods, and sequences. Your files can be located here or you can locate these files anywhere that is accessible to these programs.



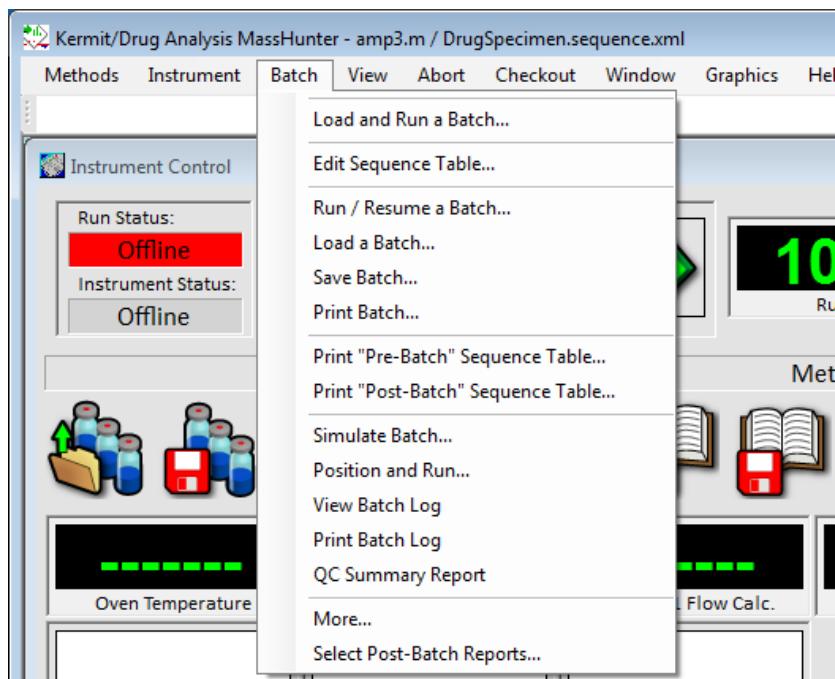
- **The data directory** contains the data from each batch run, stored in an individual directory.
- **The methods directory** contains all of your methods. Each method is stored in its own (.m) subdirectory, which contains all of the files for the method, such as compound lists, calibration tables, reports, etc.
- **The sequence directory** contains all of your sequence files. (The MassHunter GCMS Acquisition format sequence files use the (.sequence.xml) extension.)

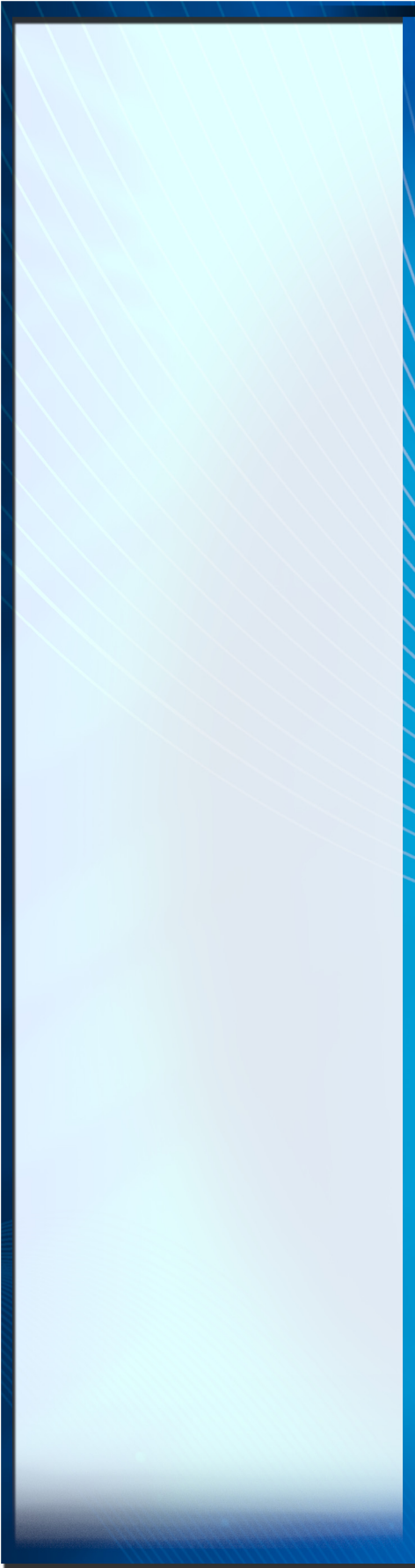
Review the Batch Menu

The few menu items that are unique to MassHunter GCMS Acquisition when it is configured in the Drug Analysis Workflow mode are described in this section.

When MassHunter GCMS Acquisition is configured in the Drug Analysis Workflow mode, the **Batch** menu contains the options associated with loading, editing, and running batches and sequences, plus access to Drug Analysis QC reports.

When MassHunter GCMS Acquisition is configured in the Enhanced Workflow mode, this menu is labeled **Sequence**, but contains similar options.





2

About Intelligent Sequencing

What is Intelligent Sequencing? 16

How does it work? 16

Where are limits and parameters defined? 16

What actions can be taken by Intelligent Sequencing? 18

What are some examples of how Intelligent Sequencing can be used? 19

How are Blanks handled? 20

How are Negatives and Controls handled? 24

How are Specimens handled? 30

Maximum # of Retries 37

Review the Intelligent Sequencing Feature

Learn How it Works

Review Examples

Understand How Blanks, Negatives, Controls, and Specimens are Handled



Agilent Technologies

What is Intelligent Sequencing?

How does it work?

Where are limits and parameters defined?

Definition - Intelligent Sequencing capabilities provided in the Drug Analysis Workflow mode of MassHunter GCMS Acquisition, automatically adjust the sequence it is running based on the quantitative and qualitative results of each analyzed sample.

Intelligent Sequencing will analyze each sample as it is processed, then, if that analysis does not meet the criteria set for your analysis, the system will perform the corrective actions defined in Intelligent Sequencing. For example it could, reinject the sample, inject a blank, skip to the next batch, or pause the sequence.

Generally speaking, when processing a sequence using the Intelligent Sequence option the following steps occur.

- 1 The sequence table identifies the samples to be run, what *Type* of samples they are (i.e., Blank, Negative, Control, or Specimen), and what method to use to process each sample.
- 2 When the sequence begins, the data for the first sample is collected and analyzed.
- 3 The results of the analyzed sample are compared to the criteria limits you specify in your method.
- 4 **If the data are within your acceptable limits**, the next sample is processed.
- 5 **If the data are outside your acceptable limits**, the decisions specified in the Intelligent Sequencing portion of the method are used to continue or pause your batch. Using those instructions, the system may:
 - a Inject a blank before continuing with the next sample in the sequence
 - b Pause the sequence
 - c Re-inject the sample
 - d Skip to the next batch
 - e Etc...

The criteria you use, and therefore, the decisions Intelligent Sequencing makes, may be based on agency regulations, good laboratory practices, or simply the unique needs of your laboratory.

Intelligent Sequencing analyzes each of the different sample types in a different way, based on the criteria you specify in the method. When you create your method, these criteria are entered in two places.

- In the Data Analysis portion of the Method you will set the acceptance criteria limits for Internal Standards, Controls, Negatives, and Blanks. See [“Step 7: Set up outliers for intelligent sequencing decisions” on page 61](#).
- In the Data Acquisition portion of the Method you will set the Intelligent Sequencing Parameters as described in the next section.

Intelligent Sequence Parameters

Here you will define how you want the sequence to proceed if an analyzed sample (Blank, Negative, Control, or Specimen) falls outside the acceptable limits you specified.*

Details for how to complete the entries for each of the 4 sample types, highlighted in this example, are provided on the following pages:

- Blanks (See [page 21](#))
- Negatives (See [page 25](#))
- Controls (See [page 25](#))
- Specimens (See [page 31](#))

You may access this dialog box at any time by selecting **Methods/Intelligent Sequencing Parameters**.

*See “[Step 7: Set up outliers for intelligent sequencing decisions](#)” on [page 61](#) for details on how to set up your acceptable limits for these sample types.

This dialog is unique to MassHunter GCMS Acquisition configured for the Drug Analysis Workflow mode.

Intelligent Sequencing Parameters

Criteria for Blanks		Criteria for Controls	
First Blank vial	91	If ISTD or analyte criteria are not met	Reinject a Control
Number of Blanks	5	If concentration is not correct	Inject a blank before reinject
Inject a blank every	3 injection(s)	Maximum # of reinjections	3
Maximum # of reinjections	3		
If one blank is contaminated	Inject blank from next vial		
		Criteria for Specimens	
If all blanks are contaminated, batch will pause.		If ISTD or analyte criteria are not met	Reinject a specimen
		If concentration is > carryover level	Inject a blank before reinject
Criteria for Negatives		<input type="checkbox"/> Enable Small Inject of:	0.00 μ L
If ISTD criteria are not met	Reinject a negative	If chromatographic checks fail	Ignore / Continue
If analyte(s) found in negative	Inject a blank before reinject	If calibration has expired	Ignore / Continue
Maximum # of reinjections	2	Calibration Time Limit (hours)	24
		Maximum # of reinjections	2
		If Maximum # of retries reached	Ignore / Continue

OK Cancel Help

What actions can be taken by Intelligent Sequencing?

As each sample is analyzed with Intelligent Sequencing, the sample results will be used to modify the next sequence action if some, predefined, criteria are met. The Table below shows the possible sequence modifications that can occur with Intelligent Sequencing.

- If they are within the acceptable limits, the sequence continues.
- If they are outside the acceptable limits, Intelligent Sequencing can take one of the actions shown in the table below.

Acceptable Limits are defined in the Quantitative Data Analysis method. See “[Step 7: Set up outliers for intelligent sequencing decisions](#)” on page 61 for details on setting acceptable limits.

Actions to take are defined in the Intelligent Sequencing dialog presented in the Intelligent Sequencing portion of your method. See the pages listed below for more details on completing the entries for each of these 4 sample types:

- Blanks (See [page 20](#))
- Negatives (See [page 24](#))
- Controls (See [page 24](#))
- Specimens (See [page 30](#))

Actions available based on the sample type	Sample type			
	Negative	Control	Specimen	Blank
Reinject ... Reinjects the Negative, Control, or Specimen and then tests those results before continuing.	✓	✓	✓	
Inject a blank before reinject Injects a blank before the system reinjects the specimen.	✓	✓	✓	
Ignore/Continue Allows the system to ignore the result and proceed with remainder of batch sequence.	✓	✓	✓	
Jump to Next Batch Jumps to the next line in the Sequence Table labeled with the keyword NewBatch . If there is none, the batch will pause.	✓	✓	✓	
Inject a blank before continuing Ensures that the GC/MS system is flushed with a blank before it injects the next vial.			✓	
Pause batch Pauses and waits for operator intervention either to end or continue the sequence. The injection counter for blanks is reset whenever a batch is paused.	✓	✓	✓ *	✓
Inject blank from next vial Injects a blank from the next blank vial position and continues with remainder of the sequence.				✓

* When calibration for a specimen has expired, the batch can be paused.

What are some examples of how Intelligent Sequencing can be used?

Example 1: Carryover Limit Exceeded

Intelligent Sequencing allows quality guidelines to be imposed without operator input. Here are a few examples.

Instead of having an operator examine every injection as it occurs and manually injecting a blank when needed, you may set the criteria which specifies:

If a specimen is found to contain a drug above the carryover limit, automatically inject a blank before continuing with the next sample in the Sequence Table.

This can save you valuable laboratory time by:

- not having to re-extract specimens contaminated by carryover
- not having to rerun specimens contaminated by carryover
- allowing the instrument operator to perform other duties

Example 2: Drug Found in the Negative

With Intelligent Sequencing if a drug of interest is found in the negative sample, the batch will be stopped early and corrective action can be taken before the bad negative invades the entire batch.

Consider the case when a negative is found to contain the drug of interest above the maximum concentration allowed for negatives. If Intelligent Sequencing is not used in this situation (and the instrument operator is not present to make the appropriate decision), then the batch will continue to run. The bad negative will invalidate the entire batch. Thus, all samples in the batch sequence will have to be reextracted and rerun. In this case you may set Intelligent Sequencing criteria to:

“Pause if the drug of interest is found in the negative.”

Important Considerations

Intelligent Sequencing decisions need to be set with care. If, for example, a negative is found to contain the drug of interest, and the decision criteria are set to “inject a blank before reinjection” and the number of reinjections is set to 9, the run may be as follows:

- original injection of the negative,
- a blank injection (because drug was found in the negative),
- a reinjection of the negative (because you set Intelligent Sequencing to re-inject),
- followed by a blank injection (if the re-injected negative still contained drug)
- and a negative rejection (if each re-injected negative contained drug)

This run would be repeated 8 more times for a total of 19 injections, wasting valuable time. Thus, it is important to set the Intelligent Sequencing decisions very carefully.

How are Blanks handled?

Definition of a Blank Sample Type

For DrugQuant, a Blank is a sample that does not contain any analytes or ISTD. A Blank is generally made up of a solvent and does not normally undergo the same extraction and preparation procedures as the other sample types. This sample type can, for example, be used in a sequence between valid specimens to evaluate carry-over contamination from a previous sample, or to flush the system before injecting another sample.

You can physically place a blank sample in any tray of the autosampler, or you can use the Intelligent Sequencing feature which injects a blank at your predefined settings.

How Intelligent Sequencing Processes Blanks

Before processing any sample, the system first checks to see if it is time to inject a blank. If the injection count indicates it is time to inject a blank, the blank is injected and processed before this sample is processed. If the count is below the blank inject value the system injects the next sample and then increments the blank injection counter.

When a Blank is injected, data are acquired by the MSD and sent to MassHunter Quantitative Analysis for analysis.

MassHunter Quantitative Analysis checks the analysis to verify the results are within your Blank sample acceptance criteria. For a Blank, this includes:

- Finding a specified target compound in a concentration above what is specified in the Carryover Amount outlier

Or

- Finding a specified ISTD in a concentration greater than what is specified in the Blank Response outlier

See [“Step 7: Set up outliers for intelligent sequencing decisions”](#) on page 61.

If the analysis shows that the results are within your specified limits:

- a The Blank cycle counter is initialized. This counter is used by Intelligent Sequencing to keep track of the number of non-blank injections allowed before automatically injecting a blank.
- b The next sample in the batch is processed.

If the analysis indicates that the results are outside your specified limits, the system processes the sample based on the parameters you set in the Intelligent Sequencing parameters dialog (shown below).

[Figure 1](#) on page 23 shows a decision flow chart for a Blank sample.

Intelligent Sequencing Parameters for Blanks

Highlighted fields are described below.

1. Where the first Blank vial is located in your autosampler tray. (Position **91** in this case.)
2. How many contiguous Blank vials you have on the tray (1-10). (This example is showing **5**. So locations 91 through 95 contain Blank vials.)
3. When you want a Blank injected. (**After every 3** injections, in this example.) See “[Example – Effects of resetting the Blank counter](#)” on page 22 for more details.)
4. The number of times the system can consecutively reinject a Blank when the blank sample analysis is out of spec (0-9). (This example shows **3**.)
5. What to do if one blank is contaminated. Options include:

Pause batch - Waits for operator intervention.*

Inject blank from next vial - Gets the next Blank, makes the injection and does the analysis once again. If there are no more Blank vials, the batch is paused and the system waits for operator intervention.

The first group box in the Intelligent Sequencing dialog is dedicated to setting the Intelligent Sequencing parameters for Blanks. You may access this dialog in MassHunter GCMS Acquisition through the **Edit Entire Method** option, or by selecting **Methods/Edit Intelligent Sequencing Parameters**.

Here you will define the Blank vial locations and capacity, the default frequency of blank injection and the actions of the sequence if the blank is contaminated.

The screenshot shows the 'Intelligent Sequencing Parameters' dialog box. The 'Criteria for Blanks' section is highlighted with a red box and numbered 1 through 5. The parameters are as follows:

Criteria for Blanks	Value
First Blank vial	91
Number of Blanks	5
Inject a blank every	3
Maximum # of reinjections	3

Additional parameters in the dialog include:

- Criteria for Controls:**
 - If ISTD or analyte criteria are not met: Reinject a Control
 - If concentration is not correct: Inject a blank before reinject
 - Maximum # of reinjections: 3
- Criteria for Specimens:**
 - If ISTD or analyte criteria are not met: Reinject a specimen
 - If concentration is > carryover level: Inject a blank before reinject
 - Enable Small Inject of: 0.00 µL
 - If chromatographic checks fail: Ignore / Continue
 - If calibration has expired: Ignore / Continue
 - Calibration Time Limit (hours): 24
 - Maximum # of reinjections: 2
- Criteria for Negatives:**
 - If ISTD criteria are not met: Reinject a negative
 - If analyte(s) found in negative: Inject a blank before reinject
 - Maximum # of reinjections: 2
- Global:**
 - If Maximum # of retries reached: Ignore / Continue

Criteria defining what constitutes a Contaminated Blank are set in the Data Analysis portion of the method. See “[Step 4: Complete the Intelligent Sequencing parameters](#)” on page 42 for details.

* The injection counter for Blanks is reset whenever a batch is paused.

Automatic reset of the Blank cycle counter

Be aware:

- The Blank cycle counter resets to zero after ANY Blank is run. Blanks may be injected at times other than those specified by the default frequency. For example, you may have Intelligent Sequencing inject a blank automatically when your criteria are not met for Negatives, Controls, and Specimens. Also, you may choose to physically place blanks in any location in the autosampler tray. When these blanks are processed, similar to when a blank is injected based on the default frequency setting, the blank cycle counter resets back to zero. See [“How Intelligent Sequencing Processes Negative and Control Samples”](#) on page 24, and [“How Intelligent Sequencing Processes Specimens”](#) on page 30 for more details.
- Paused batches ALWAYS reset the Blank counter to zero.

Example – Effects of resetting the Blank counter

If this field is set to 4, and there is a specimen with signs of **carryover in your batch**, the Blank counter resets to zero after every 4 injections, plus after the sample that shows carryover, and as shown below.

- 1 Sample (1)
- 2 Sample (2)
- 3 Sample (3)
- 4 Sample (4)
- 5 Blank (runs automatically after 4 injections)

Blank counter resets to zero

- 1 Sample (5)
- 2 Carryover Sample (6) (Causes a Blank to be run because of the criteria you set.)
- 3 Blank

Blank counter resets to zero

- 1 Sample (7)
- 2 Sample (8)
- 3 Sample (9)
- 4 Sample (10)
- 5 Blank (runs automatically after 4 injections)

This process occurs whenever a blank is run. A Blank is run at two times:

- When Intelligent Sequencing says it is time to run a Blank
- When the Sequence Table identifies the sample being run as a Blank.

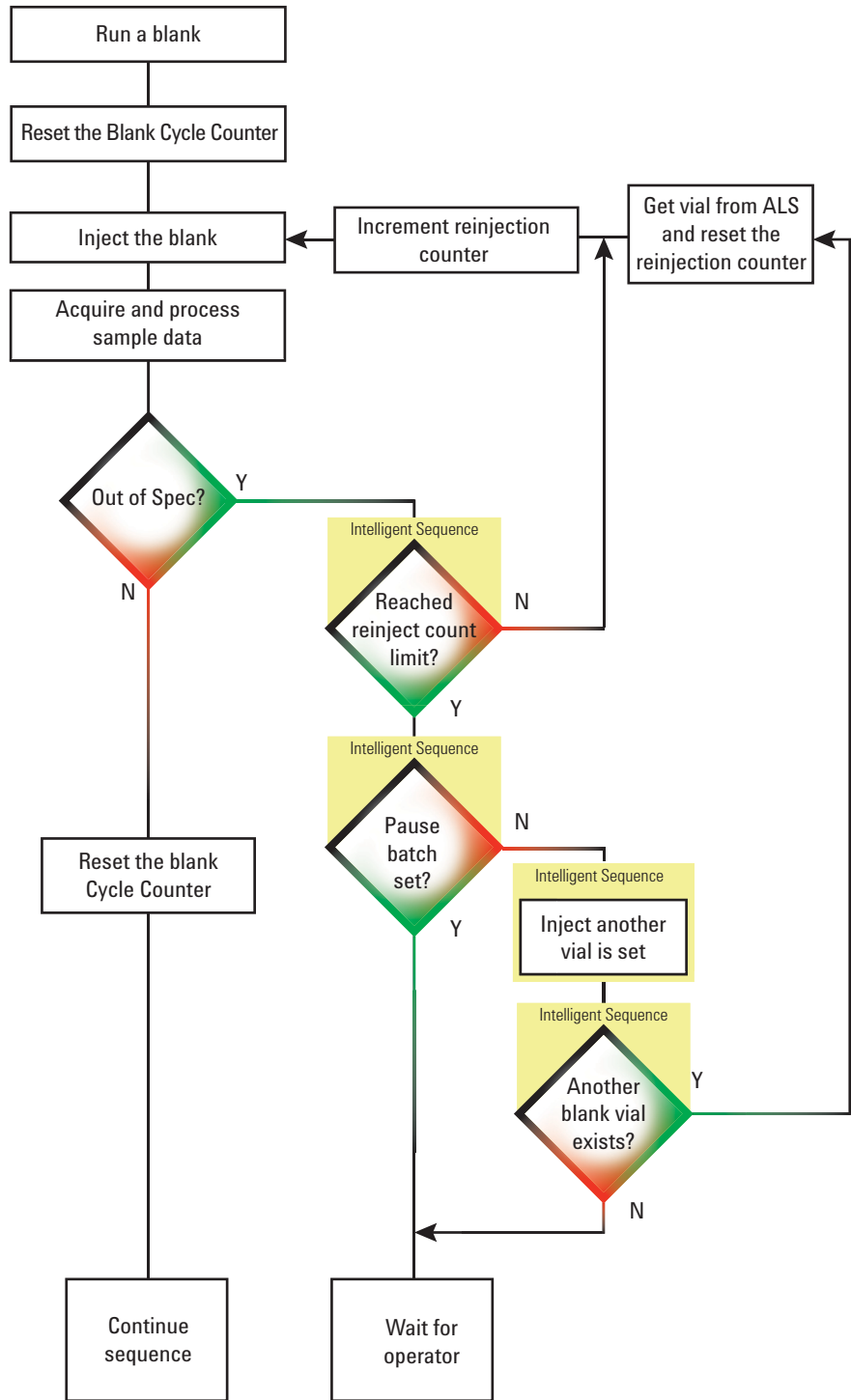


Figure 1 Blank Sample Processing Using Intelligent Sequencing

How are Negatives and Controls handled?

Definitions

How Intelligent Sequencing Processes Negative and Control Samples

A Negative Sample is a sample that contains an internal standard but no drug of interest (target compound). This sample type is generally prepared by the adding an internal standard to a clean sample matrix (urine, plasma). The negative sample goes through the same extraction and preparation procedures as all other sample types in the same batch sequence.

A Control Sample is a clean sample matrix (urine, plasma, etc.) that has been spiked with known amounts of the drug of interest and internal standard. A Control sample is a sample used in the batch to determine the integrity of the calibration and the performance of the instrument.

Before processing any sample, the system first checks to see if it is time to inject a blank. If the injection Blank cycle counter indicates it is time to inject a blank, the blank is injected and processed before this sample is processed. If the count is below the blank inject value the system injects the next sample and then increments the Blank cycle counter.

When the system processes a sample labeled as a **Negative** or a **Control**, data is acquired from the GC/MS Instrument and sent to MassHunter Quantitative Analysis for analysis. For Negative and Control samples, the processed results are checked for two things, the amount of:

- ISTD in the sample
- Drug compound ID (ion ratios) and amount in the sample

If the criteria are within your specified limits as set in your method, the system continues as usual with the next sample in the Sequence. See [“Step 7: Set up outliers for intelligent sequencing decisions” on page 61](#).

If the criteria are outside your specified limits, the system processes the sample based on the parameters you set in the Intelligent Sequencing Parameters dialog (shown below).

[Figure 2](#) on page 26 shows a flow chart of this processing.

Intelligent Sequencing Parameters for Negatives and Controls

In the **Criteria for Negatives** group box and the **Criteria for Controls** group box, you will select from dropdown lists to define what actions should be taken if the sample was shown to contain Internal Standard or Analyte(s) outside the limits you specified in your method*.

There are three areas to complete in each group. The options listed for each are shown below.

*Acceptable and unacceptable limits for Internal Standards and Analytes are set in the Data Analysis portion of the method. See “[Step 7: Set up outliers for intelligent sequencing decisions](#)” on page 61.

Criteria for Negatives	Criteria for Controls	Options include
1 If ISTD criteria are not met	1 If ISTD or analyte criteria are not met	Reinject a negative (or control). Reinjects the sample from the same vial. Pause batch. Waits for operator intervention to resolve the problem. Ignore/Continue. Ignores the result and proceeds with the remainder of the batch sequence. Jump to Next Batch. Jumps to the next line in the Sequence Table labeled with the keyword NewBatch . If there is none, the batch will pause.
2 If analytes are found in the Negative	2 If concentration is not correct	The same as those specified for the ISTD , <i>plus</i> Inject a blank before reinject. A blank is injected to remove contamination and then the sample is injected again and analyzed.
3 Maximum # of reinjections	3 Maximum # of reinjections	Here you will identify the maximum number of reinjections to make from this vial. That is, how many times you would like to re-inject and retest this sample. The range is 0 – 9. If the maximum number is reached, Intelligent Sequencing will continue based on the criteria you specify in the last field in this dialog box, labeled If Maximum # of retries reached .

This process occurs whenever a sample labeled as a negative or control is run.

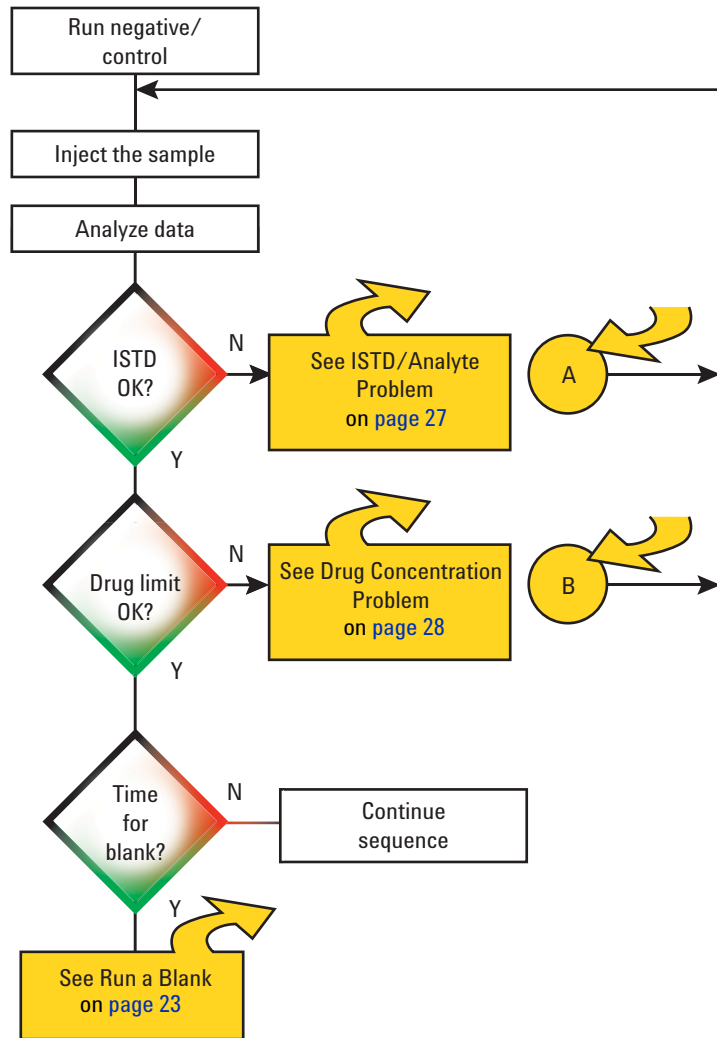


Figure 2 Negative and Control Sample Processing Using Intelligent Sequencing

This occurs whenever the internal standard is found to be outside your acceptable limits in a negative or control sample. The beginning of this process is shown in [Figure 2](#) on page 26.

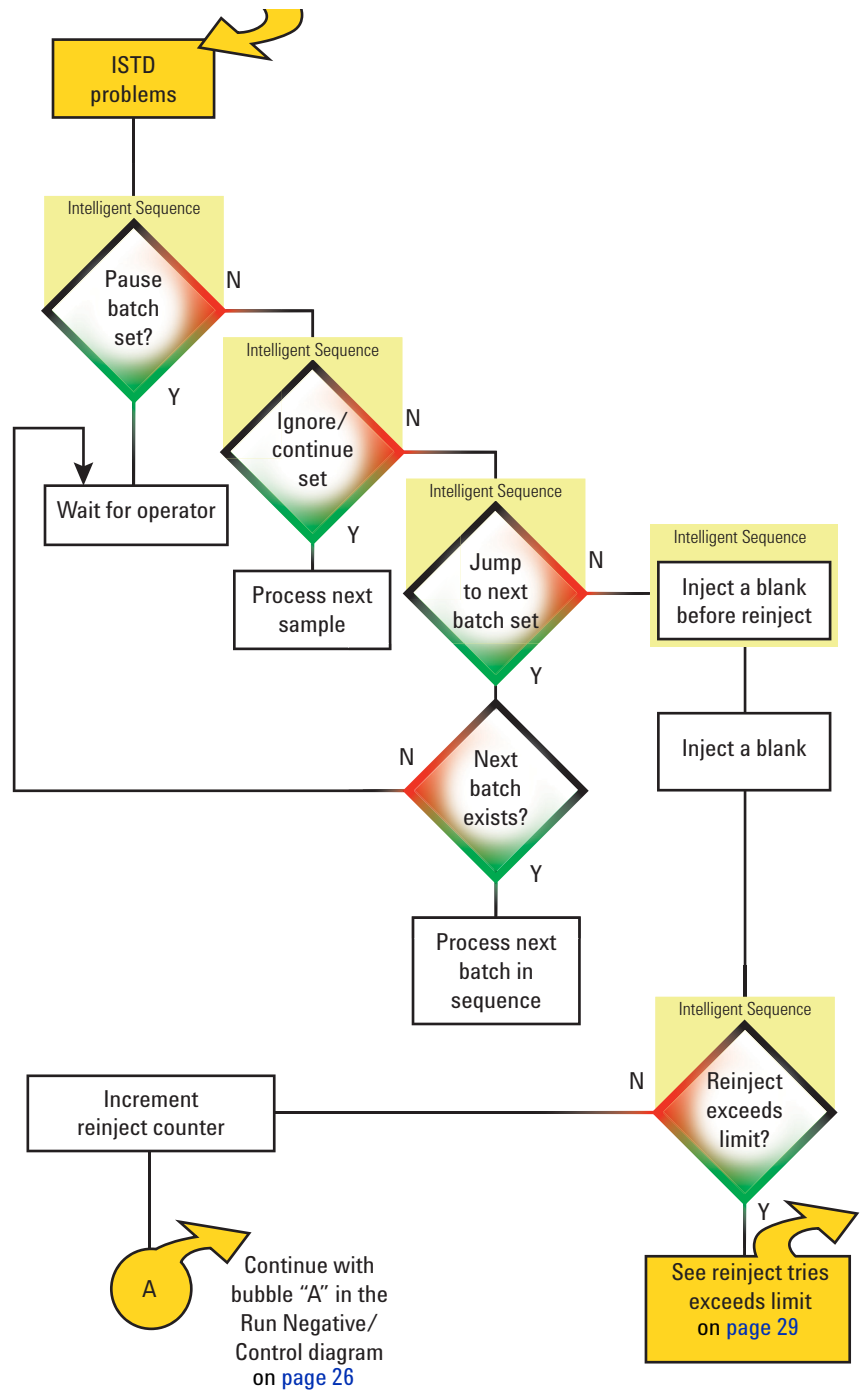


Figure 3 ISTD/analyte processing using Intelligent Sequencing

This occurs whenever the drug concentration is found to be outside your acceptable limits in a negative or control sample. The beginning of this process is shown in Figure 2 on page 26.

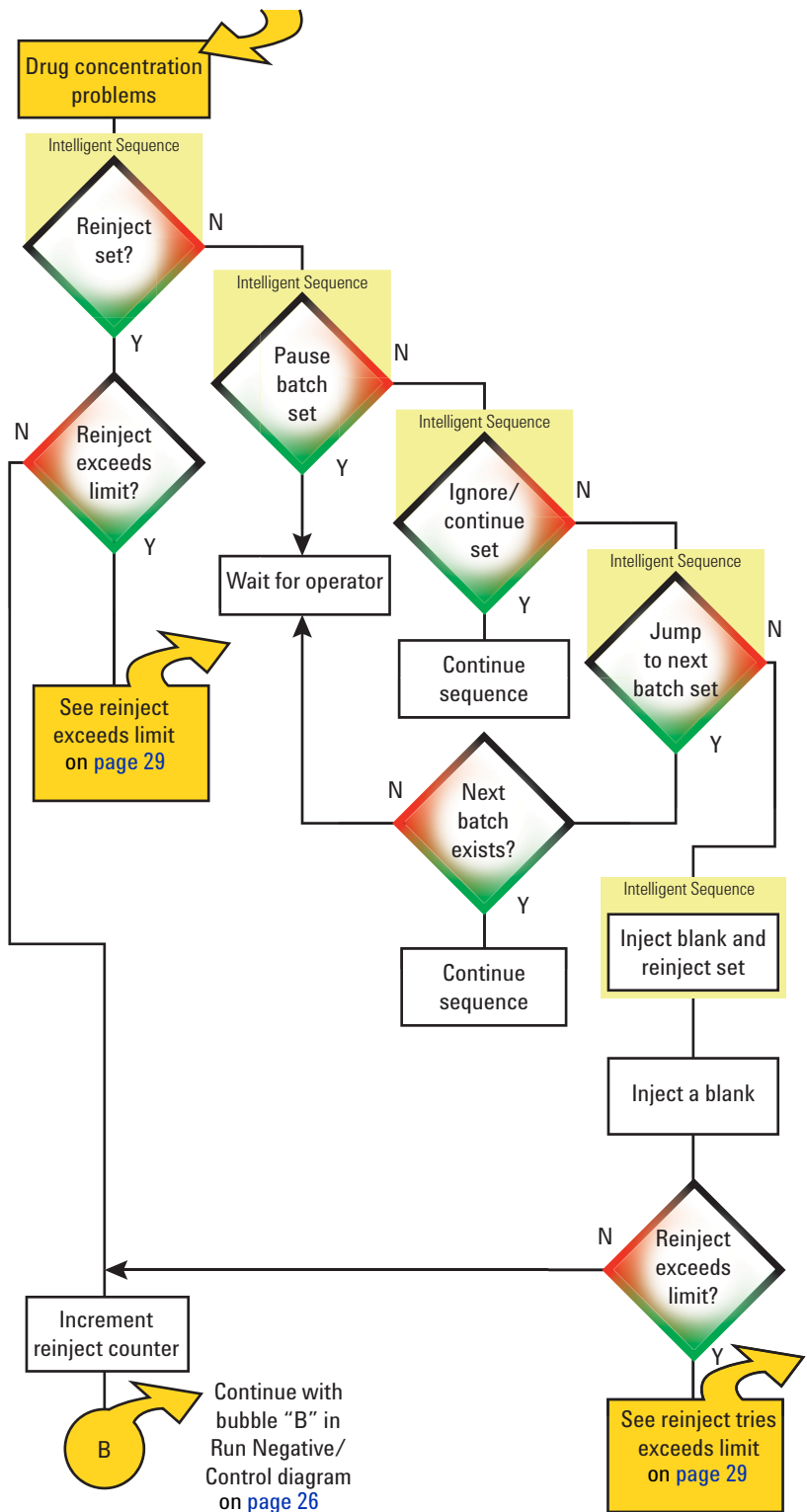


Figure 4 Drug concentration problem processing using Intelligent Sequencing

This process occurs whenever Intelligent Sequencing has specified to reinject a sample but there are no more reinjections available to process. This can happen when, for example, there are no more blank vials available for processing.

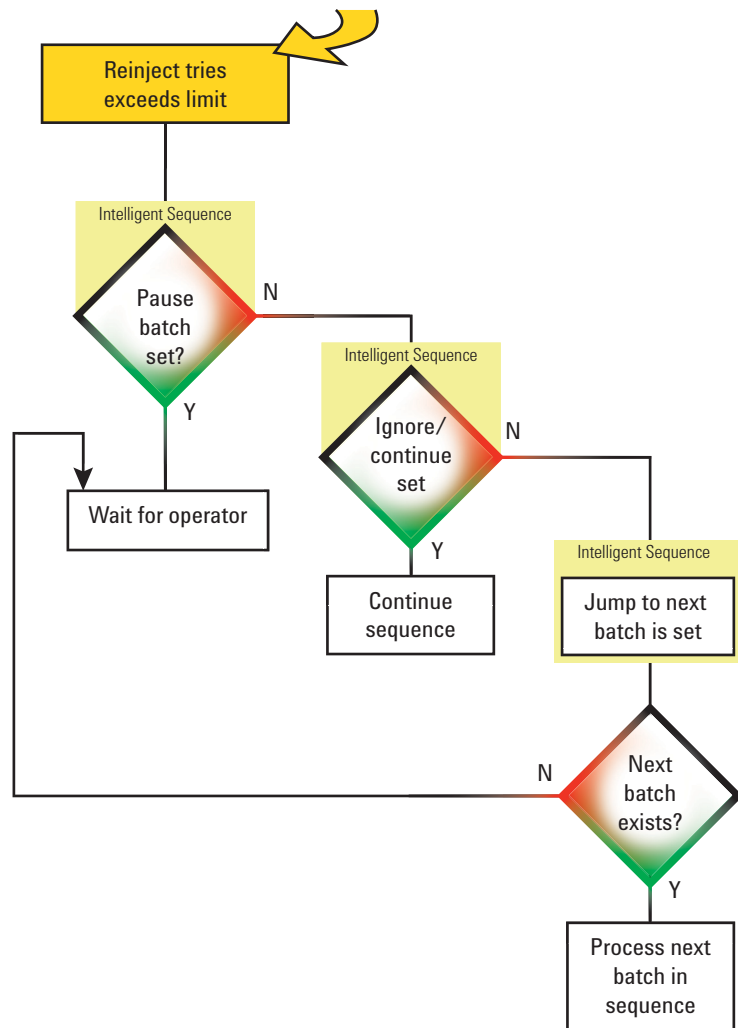


Figure 5 Reinject problem processing using Intelligent Sequencing

How are Specimens handled?

Definition of a Specimen Sample Type

How Intelligent Sequencing Processes Specimens

A Specimen sample type is a sample to be analyzed for the drug of interest.

An Internal Standard is generally spiked into the specimen and then the specimen undergoes the same extraction and preparation procedures as the other samples in the batch.

Identification of the drug of interest and the internal standard is based upon ion ratios and retention times being within your specified limits. If identification criteria are met, then the concentration of the drug of interest is calculated based on the internal standard recovery.

When the system processes a sample labeled as a Specimen, data is acquired from the GC/MS system and sent to MassHunter Quantitative Analysis for processing. For a Specimen sample MassHunter GCMS Acquisition checks the analysis for a number of items.

Checks the analysis to see if the ISTD is within the specified tolerance and the drug criteria is met. If the ISTD is outside the limits you specified in your method, the system processes the sample as described in the Intelligent Sequencing Parameters settings. [Figure 7](#) on page 33 shows a flow chart of this process.

Checks to see if the concentration of drug is greater than the carryover level allowed. The concentration is compared to maximum allowable concentration entered for the method. If it is greater than this value, it proceeds as specified in the Intelligent Sequencing parameters. [Figure 8](#) on page 34 shows a flow chart of this process.

After checking the concentration level, the system checks the **Chromatographic parameters** in the method for minimum signal to noise, resolution, fronting, tailing, and minimum/maximum peak width. If any of these parameters are out of your specified limits the system proceeds as specified in the Intelligent Sequencing parameters. [Figure 9](#) on page 35 shows a flow chart of this process.

Then the system checks to see if **Calibration time** limit in hours has expired according to limit set in Intelligent Sequencing Criteria for Specimens. [Figure 10](#) on page 36 shows a flow chart of this process.

Finally it checks the **Maximum # of reinjections** permitted. The specimen sample can only be reinjected the specified number of times set in Intelligent Sequencing Criteria for Specimens. Once that number of reinjections is reached, the system will proceed as specified in Intelligent Sequencing parameters. [Figure 5](#) on page 29.

Intelligent Sequencing Parameters for Specimens

There are 6 areas to complete here. The options for sections 1, 2, 3, and 4 are shown in the table below.

In the **Calibration Time Limit (hours)** box, enter the maximum calibration time limit, in hours. Calibrations that exceed this number of hours are considered expired. (-1 disables this parameter.) This example shows **24**. Enter what is appropriate for your method.

In the **Maximum # of reinjections** box, enter the number of reinjections to allow from this vial. That is, how many times you would allow the system to re-inject and retest this sample. The range is 0 to 9. This example show **2**. Yours may be different, depending on your method.

*Acceptable and unacceptable limits for internal standards and analytes are set in the Data Analysis portion of the method. See “Step 7: Set up outliers for intelligent sequencing decisions” on page 61.

In the Criteria for Specimens group box you will select from dropdown lists to define what Intelligent Sequencing should do when the results of a Specimen, do not meet your criteria.*

Options include	1 ISTD or analyte criteria are not met	2 Concentration is > carryover level	3 Chromatographic checks fail	4 Calibration has expired
Ignore/Continue Ignores the result and proceeds with the remainder of the batch sequence.	✓	✓	✓	✓
Reinject the specimen Reinjects the s from the same vial.	✓		✓	
Pause batch Waits for operator intervention to resolve the problem.				✓
Jump to Next Batch Jumps to the next line in the Sequence Table labeled with the keyword NewBatch . If there is none, the batch will pause.				✓
Inject a blank before reinject A blank is injected to remove contamination and then the sample is injected again and analyzed.		✓		
Inject a blank before continuing A blank is injected to remove contamination and then the next sample in the batch is injected and analyzed.		✓		

This process occurs each time a specimen is processed.

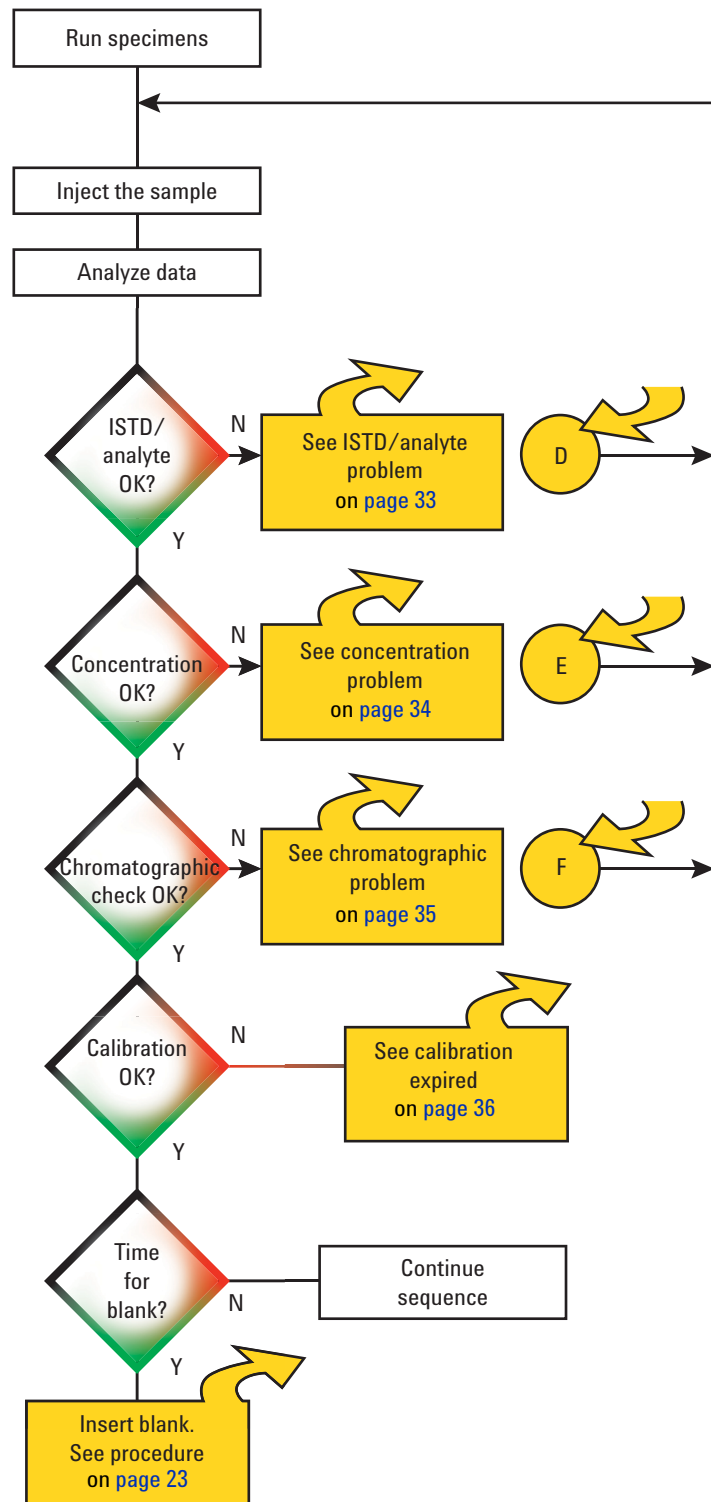


Figure 6 Specimen Sample Processing Using Intelligent Sequencing

This process occurs when an internal standard or analyte found in a specimen is outside your acceptable limits.

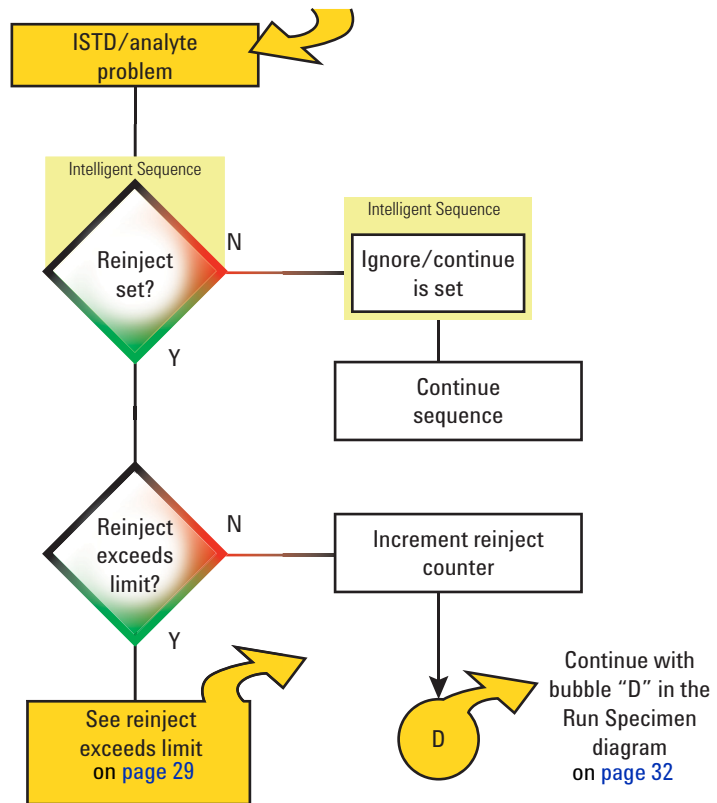


Figure 7 ISTD/analyte processing using Intelligent Sequencing

This process occurs when the concentration of analyte found in a specimen is outside the acceptable carryover concentration range.

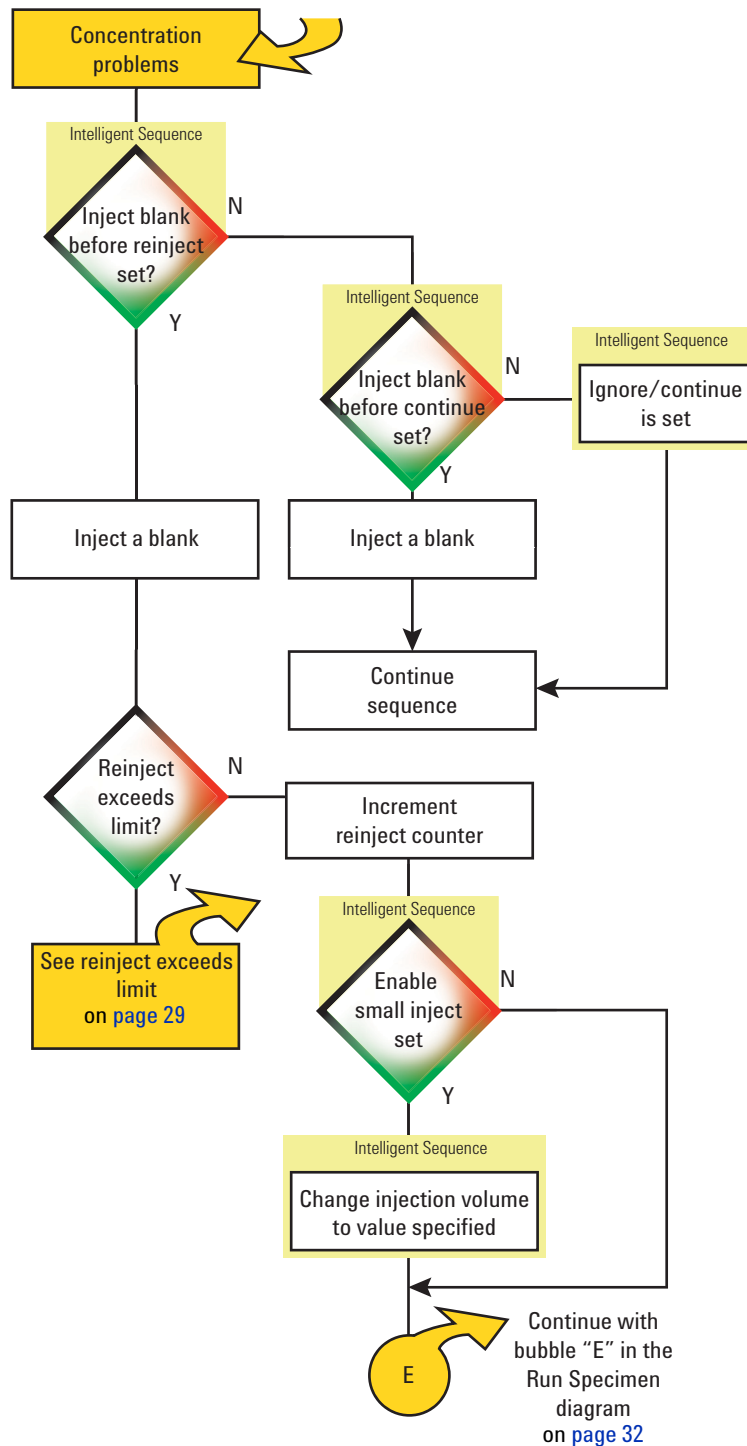


Figure 8 Drug concentration problem processing using Intelligent Sequencing

This process occurs when the signal to noise, peak resolution, peak symmetry, or peak width of a chromatogram for a specimen are outside your acceptable limits.

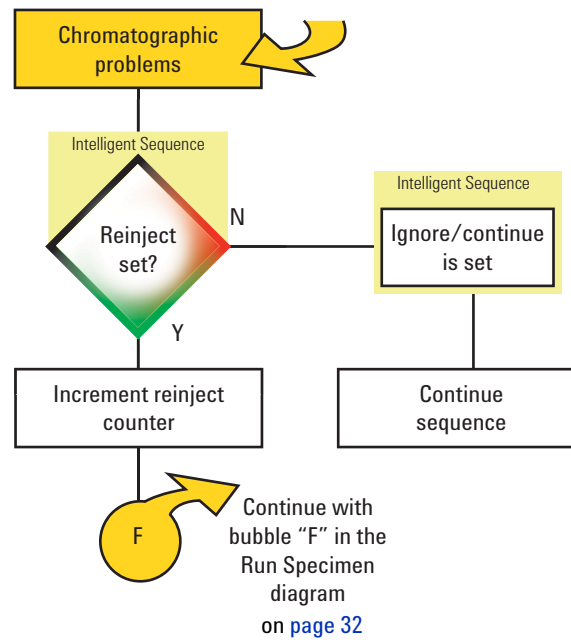


Figure 9 Chromatographic problem processing using Intelligent Sequencing

This process occurs if, when running a specimen, the system recognizes the calibration has expired.

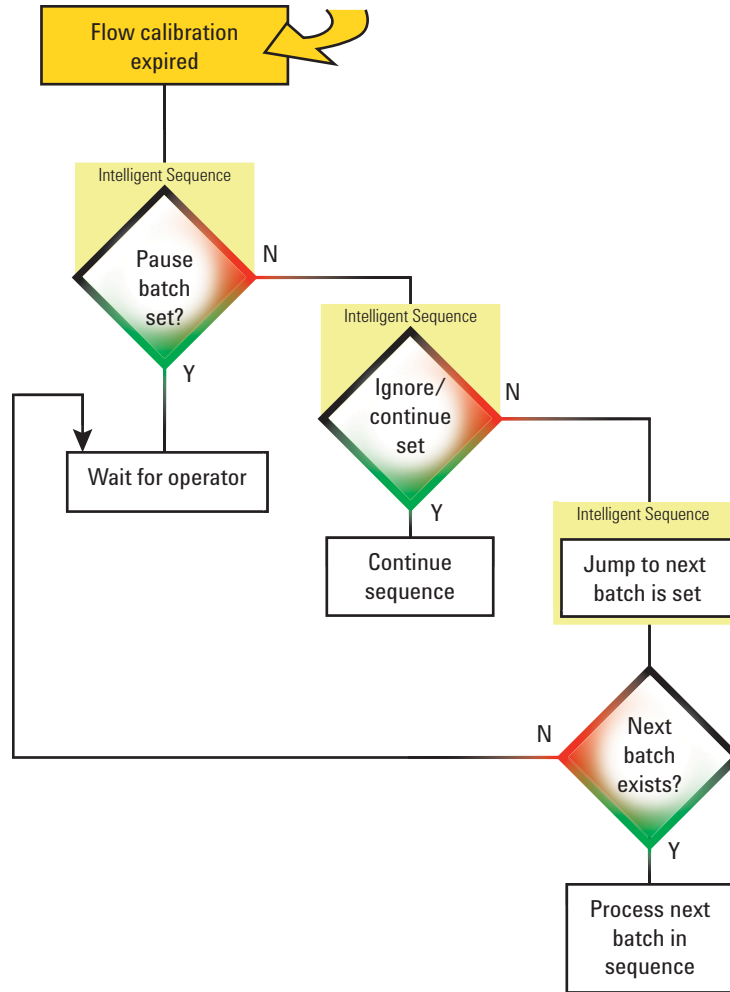


Figure 10 Calibration expired problem processing using Intelligent Sequencing

Maximum # of Retries

The last box on the Intelligent Sequencing Parameters dialog is the **Maximum # of retries reached** box.

Here you will define what to do when any of the “**Maximum # of reinjections**” options is reached for:

- Blanks
- Negatives
- Controls
- Specimens

From the drop-down list box, select:

- **Ignore/continue** – which allows the system to ignore the result and proceed with the remainder of the batch sequence.
- **Pause Batch** – which pauses the batch and waits for manual intervention either to end or continue the sequence.
- **Jump to Next Batch** – Jumps to the next batch. The next batch **MUST** have been specified by the NewBatch keyword in the Sample Log Table, otherwise the batch will pause.

See [Figure 5](#) on page 29 for a flow diagram of this process.

Intelligent Sequencing Parameters

Criteria for Blanks

First Blank vial 91

Number of Blanks 5

Inject a blank every 3 injection(s)

Maximum # of reinjections 3

If one blank is contaminated
Inject blank from next vial

If all blanks are contaminated, batch will pause.

Criteria for Negatives

If ISTD criteria are not met
Reinject a negative

If analyte(s) found in negative
Inject a blank before reinject

Maximum # of reinjections 2

Criteria for Controls

If ISTD or analyte criteria are not met
Reinject a Control

If concentration is not correct
Inject a blank before reinject

Maximum # of reinjections 3

Criteria for Specimens

If ISTD or analyte criteria are not met
Reinject a specimen

If concentration is > carryover level
Inject a blank before reinject

Enable Small Inject of: 0.00 µL

If chromatographic checks fail
Ignore / Continue

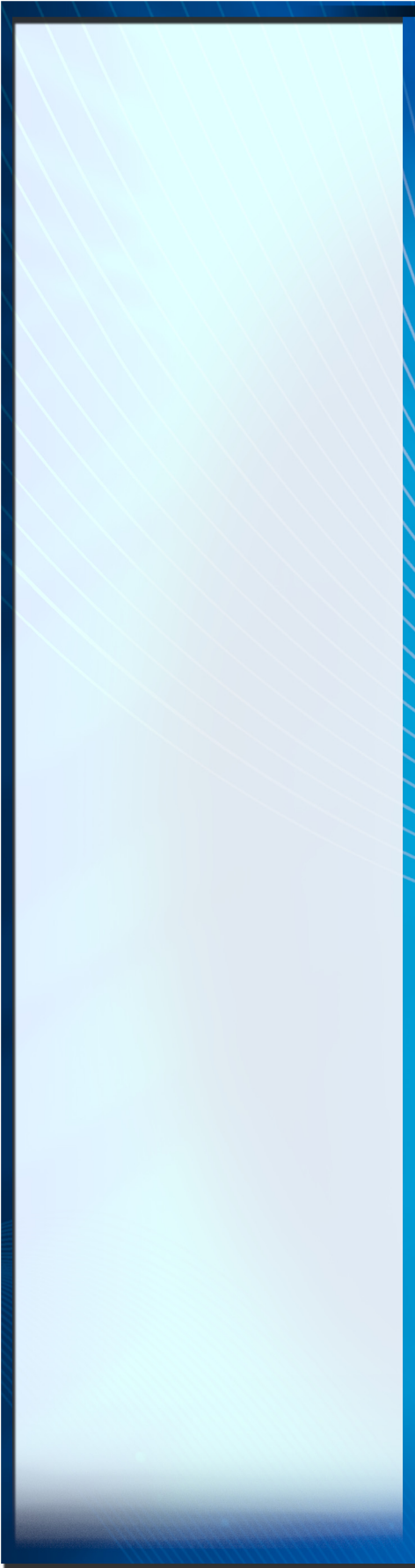
If calibration has expired
Ignore / Continue

Calibration Time Limit (hours) 24

Maximum # of reinjections 2

If Maximum # of retries reached Ignore / Continue

OK Cancel Help



3

Create the Data Acquisition Method

- Step 1: Load a data acquisition method 40
- Step 2: Select the parts of the method to edit 40
- Step 3: Describe the method and where it is saved 41
- Step 4: Complete the Intelligent Sequencing parameters 42
- Step 5: Complete the Standard Instrument Acquisition dialogs 43
- Step 6: Create a SIM method from a Scan method 44
- Step 7: Save the Method 45

Create the Data
Acquisition Method

Load a Method

Enter
Intelligent
Sequence
Parameters

Edit the
Acquisition
Parameters

Save the
Method



Agilent Technologies

Step 1: Load a data acquisition method

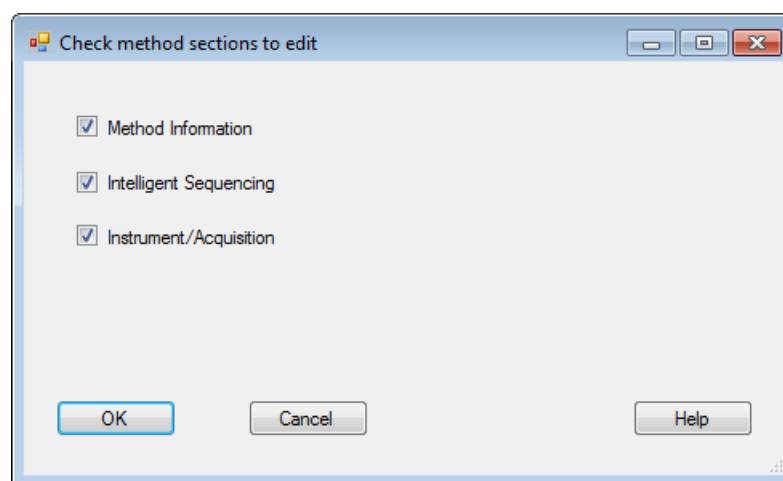
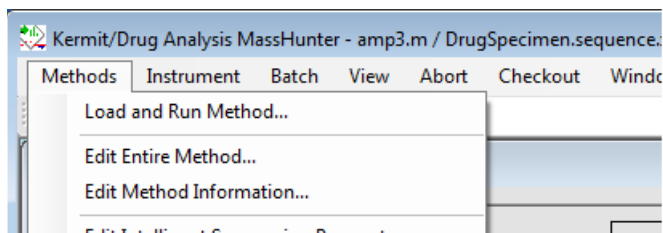
1. From the Instrument Control view menu select **Methods/ Load and Run Method...** then navigate to the optimized scan method created for your instrument, and select it.
2. Click **OK**. to close the dialog.

Step 2: Select the parts of the method to edit

1. Select **Methods/Edit Entire Method**.
2. Check each item listed.
3. Click **OK** to display the Method Information dialog.

The following procedure describes how to create a SIM data acquisition method for acquiring calibration sample data. Here we are assuming that you have previously created and optimized a scan method with good chromatographic properties that identifies the drugs target and qualifier ions and their ratios. We will use this method as a starting point and convert it to a SIM method along the way.

During this process we will cover every part of a data acquisition method.



Step 3: Describe the method and where it is saved

1. Provide a description of the method in **Method Comments**.
2. Decide whether or not to save a copy of this method in the batch data folder.
3. Select **Data Acquisition** and **Data Analysis** for the run. Although the MassHunter Quantitative Analysis method does not yet exist, you will want to run the data analysis portion of the method when it is available.

4. Click **OK** to display the Intelligent Sequencing Parameters dialog.

This is the same for all MassHunter Modes, and is described in the MassHunter Familiarization guide found in online Help. Please refer to the Familiarization guide for more details.

Method Information

Method Comments:

Analysis of amphetamine and methamphetamine and deuterated analogs as TFA derivatives from addition of MBTFA. Method is to be modified by your lab for your particular situation. Column used is DB-1701, 10m x 0.18 mm i.d., o.40 um film.]

Save Copy of Method With Data

Method Sections to Run

Pre-Run Macros/Commands

Instrument Control: Browse...

Data Analysis: Browse...

Data Acquisition

Data Analysis

Post-Run Macros/Commands

Instrument Control: Browse...

Data Analysis: Browse...

OK Cancel Help

Note – The Data Analysis method cannot be edited in the Data Acquisition program. The data analysis method can only be created or edited in the MassHunter Quantitative Analysis program.

Step 4: Complete the Intelligent Sequencing parameters

You may access this dialog box at any time by selecting **Methods/Edit Intelligent Sequencing Parameters**.

Intelligent Sequencing is unique to the MassHunter Drug Analysis Workflow mode. You will not see this dialog box unless you are running MassHunter GCMS Acquisition in the Drug Analysis Workflow mode.

Here you will define how you want the sequence to proceed if an analyzed sample (Blank, Negative, Control, or Specimen) falls outside the acceptable limits you specified.* Details for how to complete the entries for each of the 4 sample types are provided on the pages listed below:

- Blanks (See [page 20](#))
- Negatives (See [page 24](#))
- Controls (See [page 24](#))
- Specimens (See [page 30](#))

*For details on how to set up your acceptable limits for these sample types, see “[Step 7: Set up outliers for intelligent sequencing decisions](#)” on [page 61](#).

When you have completed these parameters, click **OK** to display the Inlet and Injection Parameters dialog.

Intelligent Sequencing Parameters

Criteria for Blanks

First Blank vial 91

Number of Blanks 5

Inject a blank every 3 injection(s)

Maximum # of reinjections 3

If one blank is contaminated
Inject blank from next vial

If all blanks are contaminated, batch will pause.

Criteria for Negatives

If ISTD criteria are not met
Reinject a negative

If analyte(s) found in negative
Inject a blank before reinject

Maximum # of reinjections 2

Criteria for Controls

If ISTD or analyte criteria are not met
Reinject a Control

If concentration is not correct
Inject a blank before reinject

Maximum # of reinjections 3

Criteria for Specimens

If ISTD or analyte criteria are not met
Reinject a specimen

If concentration is > carryover level
Inject a blank before reinject

Enable Small Inject of: 0.00 µL

If chromatographic checks fail
Ignore / Continue

If calibration has expired
Ignore / Continue

Calibration Time Limit (hours) 24

Maximum # of reinjections 2

If Maximum # of retries reached Ignore / Continue

OK Cancel Help

Step 5: Complete the Standard Instrument Acquisition dialogs

Edit the remaining Instrument Acquisition parameters. The remaining 5 dialogs in the Data Acquisition portion of the method are completed in the exactly the same way for all Work flow Modes (i.e., Enhanced, Drug Quant, Gasoline, etc.), and are described in detail the MassHunter Familiarization guide and in online Help. Please refer to that documentation for more details on those dialogs. Each time you complete a dialog's entries and click OK the next dialog is opened for edit.

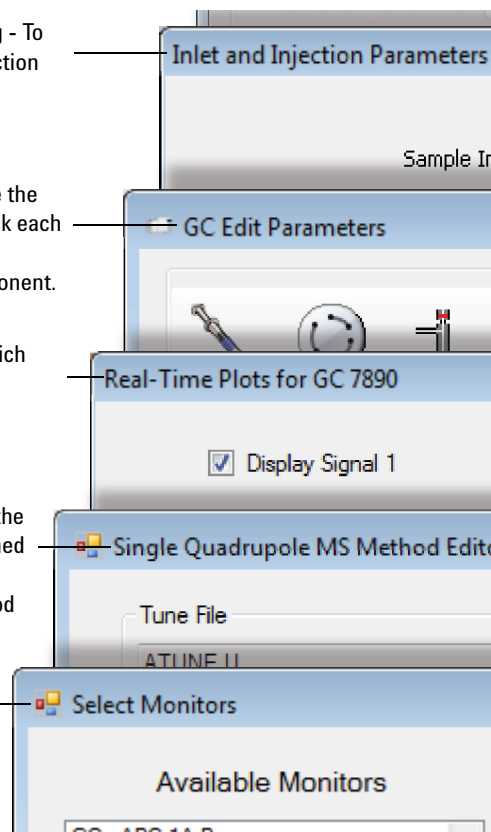
Inlet and Injection Parameters dialog - To select the sample type, inlet, and injection source.

GC Edit Parameters dialog - To define the settings for your GC. Here you will click each icon to display and complete the corresponding window for each component.

Real-Time Plots dialog - To select which signals you want displayed.

MS Method Editor dialog - To define the Tune File, SIM, Real-Time Plot, and Timed Events, settings, using the single quadrupole or triple quadrupole method editor.*

Monitors dialog - To define the MS monitors you wish to display.



*The MS Method Editor dialog is covered in more detail in “[Step 6: Create a SIM method from a Scan method](#)” on page 44.

Step 6: Create a SIM method from a Scan method

We will edit the MS Method Editor, as described here, to create an example SIM data acquisition method suitable for acquiring Amphetamine, MethAmphetamine, and their ISTDs.

Single Quadrupole MS Method Editor

Tune File: ATUNE.U

Tune Type: EI

Tune EMV: 1200

CI Gas Valve: -----

CI Flow: ----- %

MS Source: Offline 230

MS Quad: Offline 150

Run Time: 650.00 min

Solvent Delay: 1.00 min

Detector Setting: Trace Ion Detection

EM Setting: Delta EMV

Relative Voltage (V): 0.000

EM Saver: Limit Sum Limit 1e8 (Default)

Time	Start Mass	End Mass	Threshold	Scan Speed (u/s)	Frequency (scans/sec)	Cycle Time (ms)	Step Size (m/z)
1.00	50.00	550.00	500	1,562 (N=2)	2.9	342.63	0.1

Time	Group Name	Number of Ions	Total Dwell Time (ms)	Cycle Time (Hz)	Resolution	Delta EMV	Calculated EMV
1.00	Amp & methamp	8	400	2.3471	Low		1200

m/z	Dwell Time	Plot Ion
110.00	50	<input type="checkbox"/>
117.00	50	<input type="checkbox"/>
118.00	50	<input type="checkbox"/>
123.00	50	<input type="checkbox"/>
140.00	50	<input type="checkbox"/>
144.00	50	<input checked="" type="checkbox"/>
154.00	50	<input type="checkbox"/>
161.00	50	<input type="checkbox"/>

Select **SIM** as the **Acquisition Type**.

In this example SIM method, enter the ions specified here in the SIM tab as shown in the dialog:

- D6-Amphetamine (144, 123)
- Amphetamine (140, 118, 117)
- D9-MethAmphetamine (161, 123)
- MethAmphetamine (154, 118, 110)

The eight ions used in this method are specified in a single SIM time segment which has a starting time of 1 min. This means that the solvent has passed through the instrument at 1 minute. When optimizing these settings for your own method, adjust this value for the time that the solvent peak passes through your instrument.

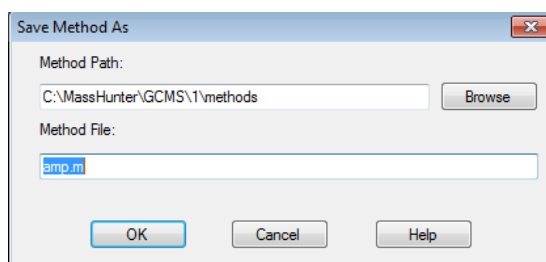
Set the Detector **EM Setting** to **Gain Factor** which is the Agilent preferred setting. The gain Factor is initially set to provide the same EMV as obtained for the most recent tune.

Step 7: Save the Method

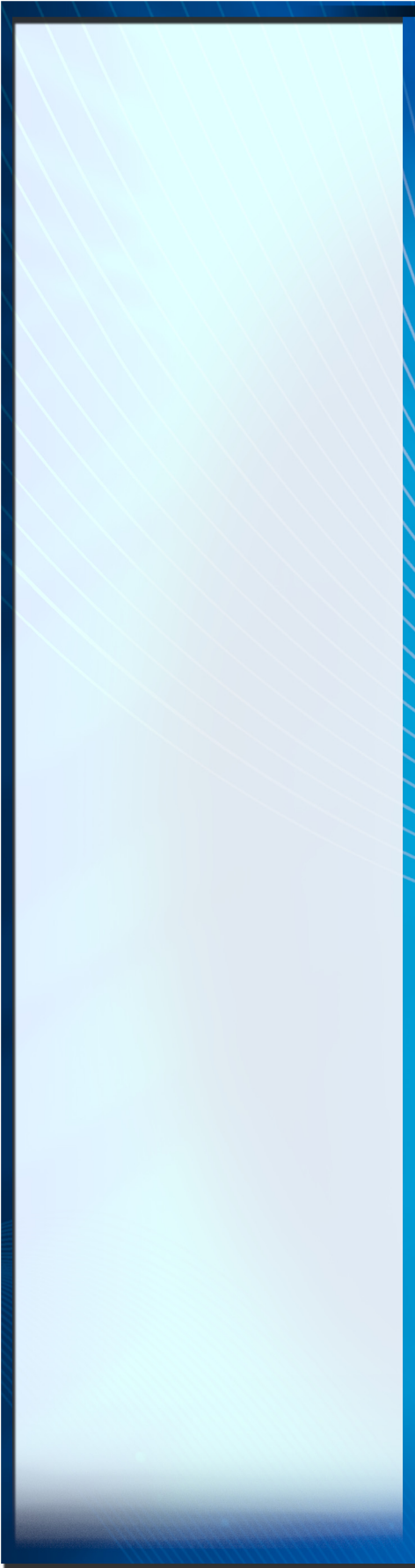
1. Enter a new name for this method.

2. Click **OK** to save the method and close the dialog.

When you have completed each of the above dialogs, Save your method.



You are now ready to continue by creating a Drug Quant method.



4

Create the Quantitative Analysis Method

Introduction	48
Step 1: Convert an MSD ChemStation method (Optional)	48
Step 2: Set up the method for Drug Quant	50
Step 3: Change each compound's Type from Scan to SIM	53
Step 4: Add a multi-level calibration to the method	54
Step 5: Add Control Samples to the Method	57
Step 6: Add ISTD Compounds to Sample as Target Compounds	59
Step 7: Set up outliers for intelligent sequencing decisions	61
Step 8: Save and Validate the Method	67

Create the Quantitative Analysis Method

Edit the Method

Set up Outliers

Save and Validate



Agilent Technologies

Introduction

Step 1: Convert an MSD ChemStation method (Optional)

1. Install the GC MSD Translator.
2. Select the starting SIM method.

In this chapter you will learn how to create a Drug Quant method that is compatible with Intelligent Sequencing.

Step 1 is an optional step. It describes how to translate an existing MSD ChemStation method into a MassHunter Quantitative Analysis method, if desired. If you are not migrating an existing MSD ChemStation method into MassHunter Quantitative Analysis, you should skip this step and begin with Step 2.

Steps 2 through 8 describe how to set up a calibration and define your acceptable limits for Blanks, Negatives, and Controls found in samples. If a sample falls outside the limits specified here, MassHunter will use the settings defined in the Intelligent Sequencing dialog to complete the sequence, as described in Chapter 2.

Also included in this chapter are details for setting up control samples used for monitoring the calibration state.

For those users interested in migrating from MSD ChemStation to MassHunter Quantitative Analysis, this step describes how to convert an existing MSD ChemStation method to a MassHunter Quantitative Analysis method using the **GC MSD Translator** tool.

If you do not want to migrate an existing MSD ChemStation method into the MassHunter Quantitative Analysis program, skip this step and create the Quant method manually by closely following the examples shown here, starting with [“Step 2: Set up the method for Drug Quant”](#) on page 50.

If the **GC MSD Translator** tool is not installed, install it using the **setup.exe** program found in the **GCMS Translator** directory located on the **Agilent MassHunter Workstation GC/MS Supplemental Software** disk.

There are 5 SIM methods supplied with your MSD ChemStation. These methods, located in the **MSDchem\msdemo\drugdemo** directory, are for:

- Amphetamines
- Cocaine
- Opiates
- PCP
- THC

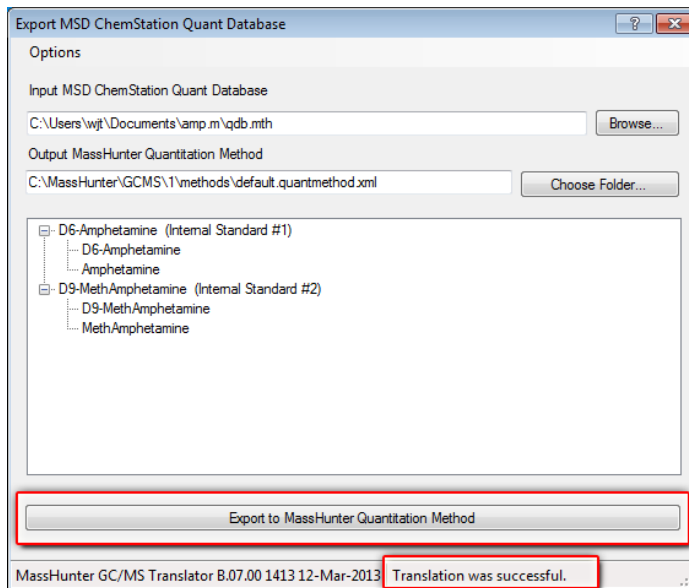
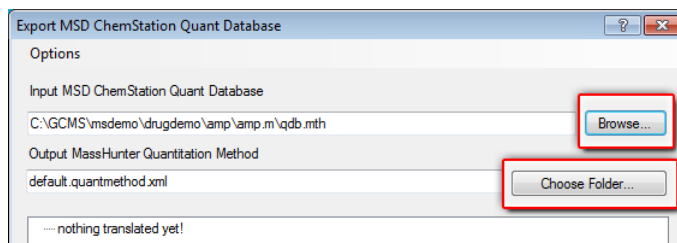
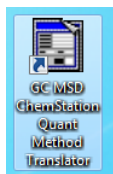
These methods use the MSD ChemStation RTE integrator in Data Analysis with customized integration parameters specific to the drugs addressed. The RTE integrator is recommended for use with MSD data when using the MSD ChemStation.

The quantitation database in these demo drug methods consists of data from a single cutoff calibration sample. The cutoff calibration sample defines the minimum concentration considered to be a positive result in a specimen sample.

The MSD ChemStation Quant Database we will use in our example is included in the **AMP.M** method, for drugs of abuse. This is a good starting point for creating a quantitative method for your analysis.

To use these methods after you convert them, you will need to edit the method parameters to suit your particular analysis requirements. You may also need to enter the calibration curve for each drug of abuse compound.

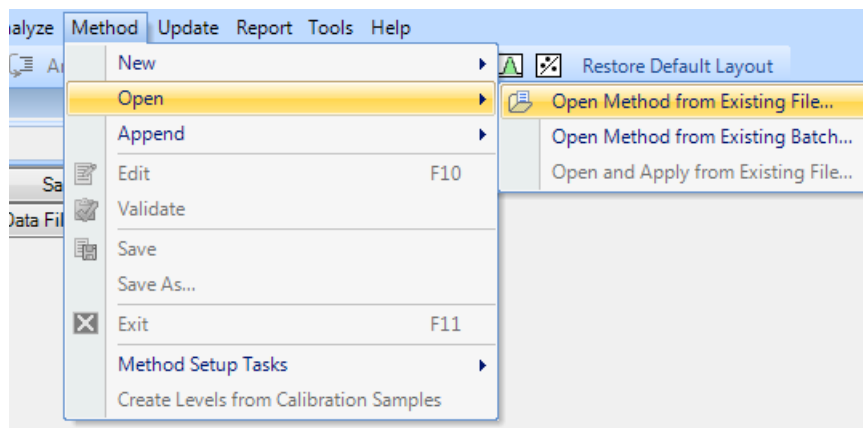
3. From the **GCMS\msdemo\drug-demo\amp** directory where your MSD ChemStation is installed, select the AMP.M method directory and copy it to a suitable location on the computer containing the **GC MSD Translator** tool.
4. Click the **GC MSD ChemStation Quant Method Translator** tool icon to start the program.
5. Click **Browse** and select the **qdb.mth** file from the MSD ChemStation method folder (i.e., AMP.M).
6. Click **Choose Folder...** and select a path for the translated file.
7. Click **Export to MassHunter Quantitation Method**.
8. If necessary, copy this new method (**default.quant-method.xml**) to the computer where the MassHunter Quantitative Analysis program is located.



Notice that the bottom right corner of the dialog now indicates **Translation was successful**. The data structure now shows the Amphetamine drugs grouped under its referenced ISTD and MethAmphetamine grouped under its ISTD.

Step 2: Set up the method for Drug Quant

1. In MassHunter Quantitative Analysis, select **Method/ Open Method from Existing File**.



If you did not convert a method, as described in the previous section, see the online Help and the Familiarization manual (accessible from online Help) for details on creating a quant method manually, then closely follow the examples in this chapter to setup the method for Drug Quant using Intelligent Sequencing.

2. Select the **default.quant-method.xml** file that you converted in the last section, or the quant method you created manually, and click **OK** to close the dialog and enter the Quant Method Editor with this method loaded.

On the next few pages, we describe how to add several levels to these calibrations so that the Target drugs will have a 4 point linear calibration curve.

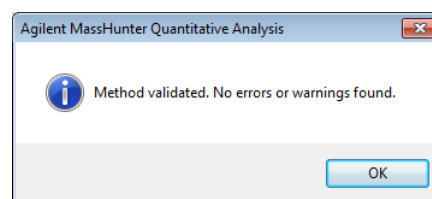
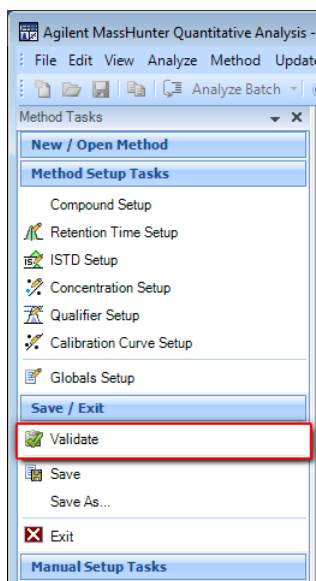
The screenshot shows the 'Method Table' window with a 'Sample' section. It contains three quantifiers, each with its own set of qualifiers and a calibration table.

Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time
Quantifier					
D6-Amphetamine		Scan	1		ISTD
Qualifier					
MZ	Rel. Resp.	Uncertainty			
123.0	65.4	20.0			
Calibration					
Level	Conc.	Response			
CU	250.0000	4995			
Quantifier					
Amphetamine		Scan	1		Target
Qualifier					
MZ	Rel. Resp.	Uncertainty			
117.0	11.7	20.0			
118.0	73.6	20.0			
Calibration					
Level	Conc.	Response			
CU	500.0000	11932			
Quantifier					
D9-MethAmphet...		Scan	1		ISTD

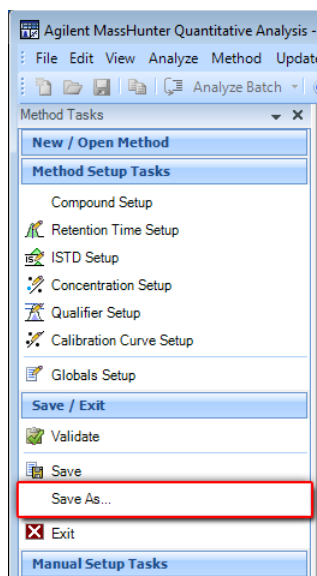
Here notice the sample structure.

- The first element is the ISTD quantifier **D6-Amphetamine**, its qualifier, and a single CU level concentration.
- Next is the **Amphetamine** Target quantifier, its qualifiers, and a single CU level concentration.

3. In the **Save/Exit** section of Method Tasks area click **Validate**. A dialog displays indicating your method is valid for MassHunter GCMS Acquisition.



4. In the **Save/Exit** section, click **Save As** and save this method as **Amp.quantmethod.xml**.

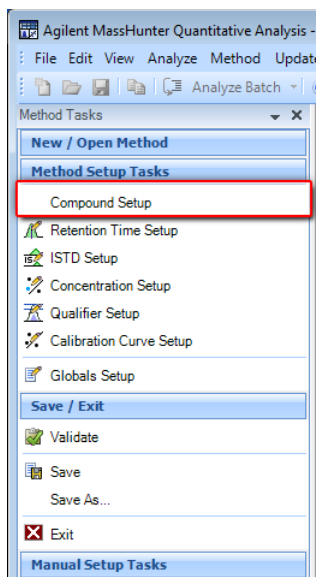


Step 3: Change each compound's Type from Scan to SIM

Skip this step if you did not convert an MSD ChemStation method.

The method translator program does not register the SIM data type automatically. To resolve this:

1. With your method loaded, in the **Method Setup Tasks** section, click **Compound Setup**.



2. In the first line of the **Scan** column select **SIM**.

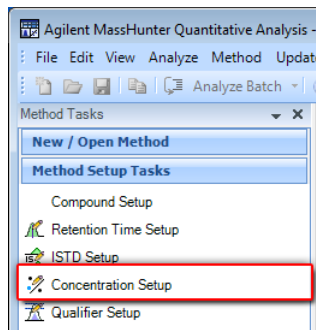
Sample					
Name	Data File	Type	Level	Acq. Metho	
Quantifier					
Name	TS	Scan	Type	MZ	
D6-Amphetamine	1	SIM	ISTD		
Amphetamine	1	Scan	Target		
D9-MethAmphet...	1	Scan	ISTD		
MethAmphetami...	1	Scan	Target		

3. Right-click on this SIM entry and select **Fill Down** from the context menu. All scans types are now correctly identified as **SIM**.

Step 4: Add a multi-level calibration to the method

1. With your method loaded, in the Method Setup Tasks section, click **Concentration Setup**.

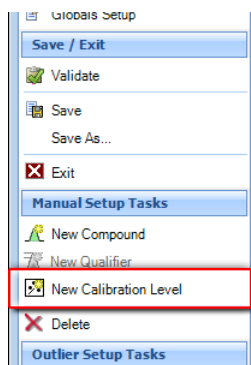
Our translated method contains a single calibration level, the CU level. In this step **we are adding 3 more calibration points** to this single point calibration to the target compounds. Later we will add control samples to this method.



2. Select the **CU Level** line in the **Amphetamine Calibration** table.

Quantifier			
Name	TS	Scan	Type
Amphetamine	1	Scan	Target
Qualifier			
MZ	Rel. Resp.	Uncertainty	
117.0	11.7	20.0	
118.0	73.6	20.0	
Calibration			
Level	Conc.	Response	
CU	500.0000	11932	

3. In the Manual Setup Tasks section, click the **New Calibration Level** icon **3 times** to add three lines to the Calibration table.



4. Enter your calibration levels.

For our example, shown below, we entered the level labels and concentration amount in ng/mL in the **Level** and **Conc** columns. Note that we are also changing the concentration of the CU Level to 250 ng/mL.

Quantifier				
Name	TS	Scan	Type	Units
Amphetamine	1	SIM	Target	NG/ML
Calibration				
Level	Conc.	Response	Enable	
L100	100.0000		<input checked="" type="checkbox"/>	
L400	400.0000		<input checked="" type="checkbox"/>	
L900	900.0000		<input checked="" type="checkbox"/>	
CU	250.0000	11932	<input checked="" type="checkbox"/>	

The **Response** column for all **Levels** will be updated later using calibration samples run from a sequence. For now we can use a linear response proportional to that of the **CU** level (11932/500 counts/conc).

5. Enter the Response for these levels.

Quantifier				
Name	TS	Scan	Type	Units
Amphetamine	1	SIM	Target	NG/ML
Calibration				
Level	Conc.	Response	Enable	
L100	100.0000	2386	<input checked="" type="checkbox"/>	
L400	400.0000	9545	<input checked="" type="checkbox"/>	
L900	900.0000	21478	<input checked="" type="checkbox"/>	
CU	250.0000	5966	<input checked="" type="checkbox"/>	

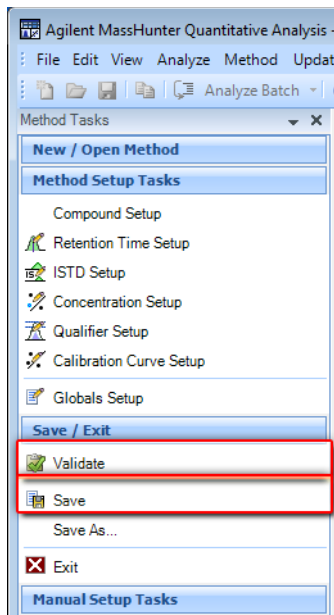
6. Use the above process to add these same levels to the Meth-Amphetamine Target quantifier (not shown).

7. Add these same levels to both ISTDs.

For each Level use an ISTD concentration of 250 ng/mL. (Only D6-Amphetamine ISTD is shown here.)

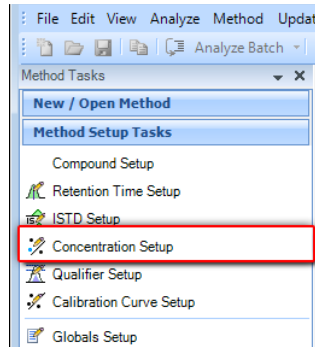
Quantifier				
Name	TS	Scan	Type	Units
D6-Amphetamine	1	SIM	ISTD	NG/ML
Calibration				
Level	Conc.	Response	Enable	
L900	250.0000	4995	<input checked="" type="checkbox"/>	
L400	250.0000	4995	<input checked="" type="checkbox"/>	
L100	250.0000	4995	<input checked="" type="checkbox"/>	
CU	250.0000	4995	<input checked="" type="checkbox"/>	

8. Validate and save the method.



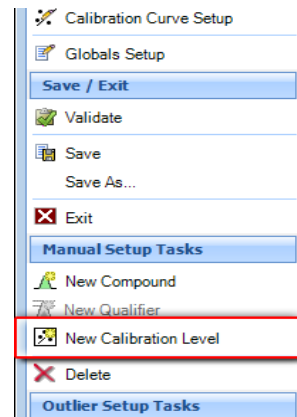
Step 5: Add Control Samples to the Method

1. With your method loaded, in the **Method Setup Tasks** section, click **Concentration Setup**.



2. Select the last line in the Calibration Table for D6-Amphetamine.

3. In the **Manual Setup Tasks** section, click **New Calibration Level** twice to add two data entry lines for new levels.



4. Fill in these levels as shown here with new controls labeled **QCLow** and **QCMed**. Since this is an ISTD the concentration will be 250 ng/mL for both.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time

Quantifier					
Name	TS	Scan	Type	Units	
D6-Amphetamine	1	SIM	ISTD	NG/ML	

Calibration				
Level	Conc.	Response	Enable	
QCLow	250.0000	4995	<input checked="" type="checkbox"/>	
QCMed	250.0000	4995	<input checked="" type="checkbox"/>	
CU	250.0000	4995	<input checked="" type="checkbox"/>	
L100	250.0000	4995	<input checked="" type="checkbox"/>	
L400	250.0000	4995	<input checked="" type="checkbox"/>	
L900	250.0000	4995	<input checked="" type="checkbox"/>	

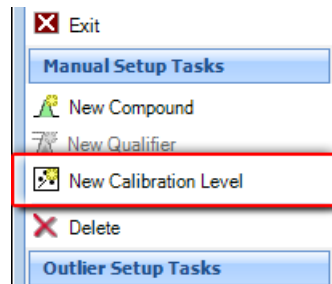
5. Use the above procedure to add these controls to the other ISTD.

6. Select the last line in the Calibration Table for Amphetamine.

7. In the **Manual Setup Tasks** section, click **New Calibration Level** twice to add two data entry lines for new levels.

8. Fill in these levels as shown here with new controls labeled **QCLow** and **QCMed**.

9. Using a procedure similar to the preceding steps, add these two controls to the other target compounds.



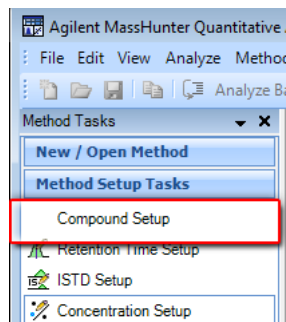
Level	Conc.	Response	Enable
QCLow	150.0000	3579	<input checked="" type="checkbox"/>
QCMed	313.0000	7468	<input checked="" type="checkbox"/>
CU	250.0000	5966	<input checked="" type="checkbox"/>
L100	100.0000	2386	<input checked="" type="checkbox"/>
L400	400.0000	9545	<input checked="" type="checkbox"/>
L900	900.0000	21478	<input checked="" type="checkbox"/>

The concentrations here are set for a regulation requiring a control set at 40% below and 25% above the cutoff (**CU**) level.

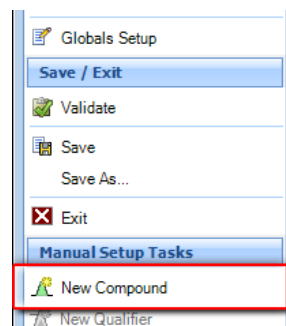
Step 6: Add ISTD Compounds to Sample as Target Compounds

1. With your method loaded, in the **Method Setup Tasks** section, click **Compound Setup**.

MassHunter GCMS Acquisition cannot process a carryover ISTD response in a Blank sample run as it does with a carryover Target compound response in a Blank. Therefore, it is necessary to set up additional compounds with the same ion identity as the ISTD we want to monitor for carryover when running a Blank sample but identify them as Target compounds.



2. In the **Manual Setup Tasks** section, click **New Compound** twice to add two lines to the Quantifier table.



3. Enter the compound Name, Scan, Type, MZ and RT, Criteria as shown here for D6Amp and D9MethAmp.

For our example, we are adding D6-Amphetamine and D9-MethAmphetamine to our method as target compounds under a different compound name.

Name	Data File	Type	Level	Acq. Method File	A		
Quantifier							
Name	TS	Scan	Type	MZ	RT	Ion Polarity	Criteria
D6-Amphetamine	1	SIM	ISTD	144.0	1.799	Positive	Close RT with Qu
Amphetamine	1	SIM	Target	140.0	1.807	Positive	Close RT with Qu
D9-MethAmphetamine	1	SIM	ISTD	161.0	1.998	Positive	Close RT with Qu
MethAmphetamine	1	SIM	Target	154.0	2.021	Positive	Close RT with Qu
D6Amp	1	SIM	Target	144.0	1.799	Positive	Close RT with Qu
D9MethAmp	1	SIM	Target	161.0	1.998	Positive	Close RT with Qu

Note that other than the name and Type, they are identical to the two ISTDs.

4. Create the Quantitative Analysis Method Step 6: Add ISTD Compounds to Sample as Target Compounds

- In the **Method Setup Tasks** section, click **Qualifier Setup**.
- Select the new compound D6Amp in the Quantifier table.
- In the **Manual Setup Tasks** section, click **New Qualifier** to add a Qualifier Table under the D6Amp quantifier with one qualifier entry line.
- Enter the **MZ, Rel. Resp.**, and **Uncertainty** for the **D6Amp** qualifier as shown here.
- Repeat similar to the above steps to add the **D9MethAmp** qualifier as shown here.
- Validate the method and save it.

For our example, we are copying the D6-Amphetamine qualifier entries here.

Qualifier				
Name	TS	Scan	Type	MZ
D6Amp	1	SIM	Target	144.0
Qualifier				
MZ	Rel. Resp.	Uncertainty	Area Sum	
123.0	65.4	20.0	<input type="checkbox"/>	

Qualifier				
Name	TS	Scan	Type	MZ
D9MethAmp	1	SIM	Target	161.0
Qualifier				
MZ	Rel. Resp.	Uncertainty	Area Sum	
123.0	25.3	20.0	<input type="checkbox"/>	

Step 7: Set up outliers for intelligent sequencing decisions

Set Limits for ISTDs in Blank Samples

- With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- Click **Blank Response** and enter the acceptable maximum response for each ISTD compound found in a Blank sample type. The example shows the entry of acceptable carryover responses for the ISTDs in a Blank sample.

Set the Acceptable Response Range for an ISTD in a Negative Sample

- With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- Click **ISTD Response** and enter the High and Low response for each ISTD compound found in a Negative sample type.

Here we set up the allowable specifications for our sample analysis. When a specification entered here is not met, a decision entered in Intelligent Sequencing is carried out by MassHunter GCMS Acquisition. Leave an entry blank if you do not want MassHunter GCMS Acquisition to report or act on that outlier.

As previously discussed in “[Step 6: Add ISTD Compounds to Sample as Target Compounds](#)” on page 59, MassHunter GCMS Acquisition cannot process an ISTD response for a Blank sample type directly. MassHunter GCMS Acquisition can process a Target compound response for a Blank sample type so we set up the D6Amp and D9MethAmp target compounds in the previous step for that purpose.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time
▶					
Quantifier					
Name	TS	Scan	Type	Max. Blank Resp.	
D6-Amphetamine	1	SIM	ISTD		
Amphetamine	1	SIM	Target		
D9-MethAmphetamine	1	SIM	ISTD		
MethAmphetamine	1	SIM	Target		
D6Amp	1	SIM	Target	10	
D9MethAmp	1	SIM	Target	10	

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time
▶					
Quantifier					
Name	TS	Scan	Type	ISTD Resp. Limit Low	ISTD Resp. Limit High
MethAmphetamine	1	SIM	Target		
D9-MethAmphetamine	1	SIM	ISTD	2000	4000
D9MethAmp	1	SIM	Target		
D6-Amphetamine	1	SIM	ISTD	2000	10000
D6Amp	1	SIM	Target		
Amphetamine	1	SIM	Target		

Set an Acceptable Carryover Concentration Range for Target Compounds

- a With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- b Click **Carryover Amount** and enter the High and Low amount in ng/mL for each compound found in any sample type.

Sample						
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time	

Quantifier						
Name	TS	Scan	Type	Amt. Limit Low	Amt. Limit High	
D6-Amphetamine	1	SIM	ISTD			
D6Amp	1	SIM	Target			
Amphetamine	1	SIM	Target	0		1E+04
D9-MethAmphetamine	1	SIM	ISTD			
D9MethAmp	1	SIM	Target			
MethAmphetamine	1	SIM	Target	0		1E+04

Set Concentration Limits for Target Compounds in Negative Samples

- With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- Click **Negative Concentration** and enter the acceptable response in ng/mL for each Target compound found in a Negative sample type.

Set an Acceptable Concentration for Target Compounds in Negatives

- With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- Click **Negative Concentration** and enter the acceptable response in ng/mL for each Target compound found in a Negative sample type.

This example shows the entry of acceptable carryover responses for the target compounds in a Negative sample.

Sample						
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time	Sample
▶						
Quantifier						
Name	TS	Scan	Type	Max. Blank Conc.		
D6-Amphetamine	1	SIM	ISTD			
D6Amp	1	SIM	Target			
Amphetamine	1	SIM	Target	5.0000		
D9-MethAmphetamine	1	SIM	ISTD			
D9MethAmp	1	SIM	Target			
MethAmphetamine	1	SIM	Target	5.0000		

This example shows the entry of acceptable carryover concentrations in ng/mL for the target compounds in a Negative sample.

Sample						
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time	Sample
▶						
Quantifier						
Name	TS	Scan	Type	Max. Blank Conc.		
D6-Amphetamine	1	SIM	ISTD			
D6Amp	1	SIM	Target			
Amphetamine	1	SIM	Target	5.0000		
D9-MethAmphetamine	1	SIM	ISTD			
D9MethAmp	1	SIM	Target			
MethAmphetamine	1	SIM	Target	5.0000		

Set the Concentration Linearity Limit

- a With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- b Click **Limit of Quantitation** and enter the lowest concentration in ng/mL for each Target compound found in a Negative sample type.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time
Quantifier					
Name	TS	Scan	Type	LOQ	
D6-Amphetamine	1	SIM	ISTD		
D6Amp	1	SIM	Target		
Amphetamine	1	SIM	Target	150	
D9-Methamphetamine	1	SIM	ISTD		
D9MethAmp	1	SIM	Target		
▶ Methamphetamine	1	SIM	Target	150	

Setup Minimum Signal to Noise for Specimen Samples

- a With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- b Click **Signal-to-Noise Ratio** and enter acceptable values for the target compounds.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time
Quantifier					
Name	TS	Scan	Type	Min. S/N	
D6-Amphetamine	1	SIM	ISTD		
D6Amp	1	SIM	Target		
Amphetamine	1	SIM	Target	10.00	
D9-Methamphetamine	1	SIM	ISTD		
D9MethAmp	1	SIM	Target		
▶ Methamphetamine	1	SIM	Target	10.00	

Setup Chromatographic Resolution for Specimen Samples

- With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- Click **Peak Resolution** and enter acceptable values for the target compounds.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time

Quantifier					
Name	TS	Scan	Type	Resolution Limit	
D6-Amphetamine	1	SIM	ISTD		
D6Amp	1	SIM	Target		
Amphetamine	1	SIM	Target	80.0	
D9-MethAmphetamine	1	SIM	ISTD		
D9MethAmp	1	SIM	Target		
MethAmphetamine	1	SIM	Target	80.0	

Setup Chromatographic Symmetry for Specimen Samples

- With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- Click **Peak Symmetry** and enter acceptable values for the target compounds. A value of 1.0 indicates perfect symmetry. A Limit Low less than 1 indicates allowable fronting and a Limit High greater than 1 indicates allowable tailing.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time

Quantifier						
Name	TS	Scan	Type	Symmetry Limit Low	Symmetry Limit High	
D6-Amphetamine	1	SIM	ISTD			
D6Amp	1	SIM	Target			
Amphetamine	1	SIM	Target	0.50	2.00	
D9-MethAmphetamine	1	SIM	ISTD			
D9MethAmp	1	SIM	Target			
MethAmphetamine	1	SIM	Target	0.50	2.00	

Set Allowable Peak Width for Specimen Samples

- a With your method loaded, in the **Method Setup Tasks** section, click **Outlier Setup Tasks**.
- b Click **Peak Full Width Half Maximum** and enter an acceptable peak width range in minutes for the target compounds.

Sample					
Name	Data File	Type	Level	Acq. Method File	Acq. Date-Time

Quantifier						
Name	TS	Scan	Type	FWHM Limit Low	FWHM Limit High	
D6-Amphetamine	1	SIM	ISTD			
D6Amp	1	SIM	Target			
Amphetamine	1	SIM	Target	0.050	3.000	
D9-MethAmphetamine	1	SIM	ISTD			
D9MethAmp	1	SIM	Target			
MethAmphetamine	1	SIM	Target	0.050	3.000	

Step 8: Save and Validate the Method

1. With your method loaded, in the **Method Setup Tasks** section click **Save / Exit**.
2. Click **Validate** to verify that your method is valid. If no problems exist click OK to acknowledge the verification message. If problems exist, a linked problem message displays in the **Method Error List** window located below the **Method Table**.
3. Click **Save As** and save this method in the directory you use for storing master methods used by automated sequences.

At this point our quantitation method contains all the settings required by Intelligent Sequencing when analyzing specimens for the presence of Amphetamines and MethAmphetamines. What remains is to replace the current compound responses that originated in our translated method with calibrated responses from the real analytical system.

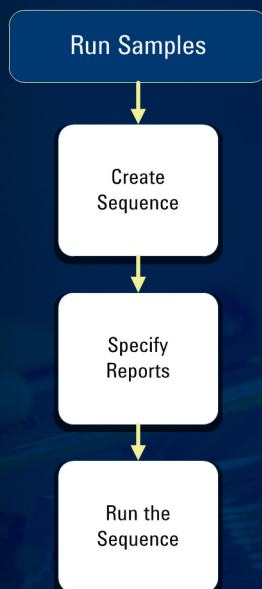
See [Chapter 6](#) for details on how to update the calibration.



5

Run Samples

Introduction 70
Step 1: Load the default Sequence 70
Step 2: Edit the Sequence Table 71
Step 3: Specify reports 73
Step 4: Save the Sequence Table 73
Step 5: Run the Sequence 74



Introduction

In MassHunter GCMS Acquisition, data acquisition is automated using the sequence table.

The Sequence Table defines the:

- Name of the sample
- Method to be used for data acquisition and analysis
- Sample Type being analyzed
- Location in which to save your results
- Optional instructions for sample processing (Keyword/Keyword string)
- ALS vial location
- Batch samples and optional name
- Reports to be run

In this chapter we explain how to set up a Sequence Table with multiple batches of samples using the keyword **NewBatch**. This keyword indicates where the MassHunter should continue processing when a *Skip to next batch* command is generated by Intelligent Sequencing.

The following example consist of HHS required Quality Control samples followed by an HHS allowable number of specimens.

Step 1: Load the default Sequence

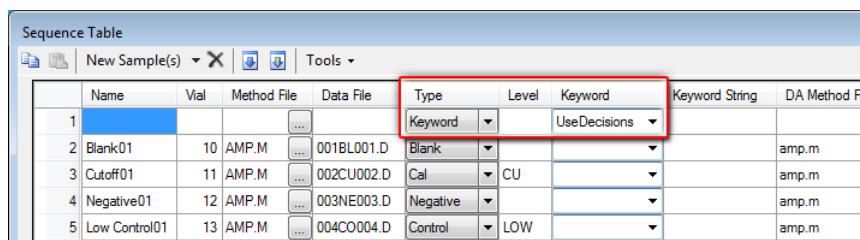
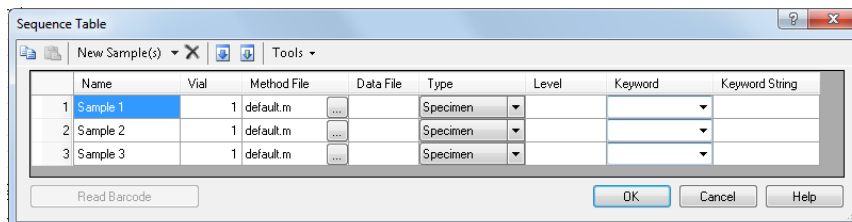
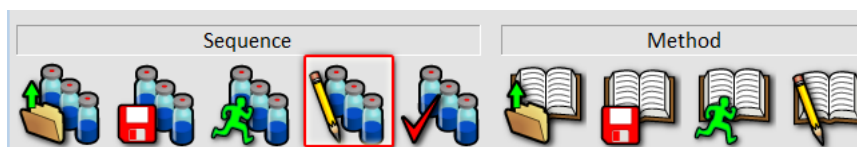
1. In the Data Acquisition Instrument Control view, click the Load Sequence icon then select the default.sequence.xml file from your instrument directory sequence folder.
2. Save this sequence as a new name in the sequence folder of the instrument directory.



If you select a different directory for storing this sequence, that location will become the default storage directory location the next time you load or save a sequence.

Step 2: Edit the Sequence Table

1. Click the Edit Sequence icon to open the Sequence Table for editing.
2. To begin, from the **Tools** menu, select **Add/Remove Columns** and add the DA Method File and DA Report Templates columns.
3. From the **Keyword** dropdown, select **Use Decisions**.



This command tells MassHunter GCMS Acquisition to send each sample's data file to MassHunter Quantitative Data Analysis for processing and wait for those results before running the next sample.

During processing, if MassHunter GCMS Acquisition finds the sample is out of spec it reviews the actions specified in **Intelligent Sequencing** and implements your previously defined decision.

In this example we are not defining the batch name (see BatchDir keyword in online help). The system names the batch directory using a unique default name consisting of date, and time, for example: 2013-06-11-1640.b. See the online help for more information.

- Start the first batch on the first line after the Keyword you just entered.

Name	Vial	Method File	Data File	Type	Level	Keyword	Keyword String	DA Method File
1				Keyword		UseDecisions		
2	Blank01	10 AMP.M	001BL001.D	Blank				amp.m
3	Cutoff01	11 AMP.M	002CU002.D	Cal	CU			amp.m
4	Negative01	12 AMP.M	003NE003.D	Negative				amp.m
5	Low Control01	13 AMP.M	004CO004.D	Control	LOW			amp.m
6	Med Control01	14 AMP.M	005CO005.D	Control	MED			amp.m

A typical sequence that includes multiple batches of specimens, starts out with a Blank sample that allows you to monitor for carryover contamination before you start to run specimens. This blank resets the Intelligent Sequencing cycle counter to zero.

For an HHS certified laboratory this is usually followed by:

- A cutoff calibration sample (**Type = CAL, Level = CU** to match the level label entered in the MassHunter Quantitative Analysis method)
- A negative sample (**Type=Negative**)
- A positive control set at 25% above the drug cutoff (**Type=Control, Level = MED**)
- A control set at or below 40% of the drug cutoff concentration (**Type=Control, Level = LOW**).

With a criteria that 10% of all samples in the batch must be quality control samples, these QC samples can be followed by up to 36 specimen samples.

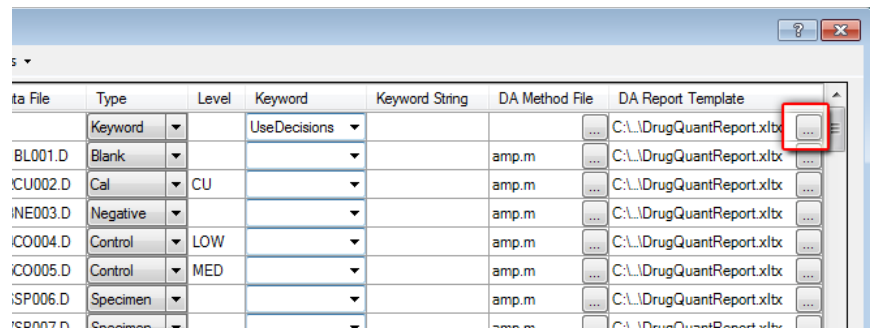
- To begin another batch in the same sequence table, set the sample type to **Keyword** and select **NewBatch** from the **Keyword** dropdown. This batch, like the first, will use the MSD ChemStation auto naming batch feature.

Name	Vial	Method File	Data File	Type	Level	Keyword	Keyword String
39	Specimen33	47 AMP.M	007SP038.D	Specimen			
40	Specimen34	48 AMP.M	007SP039.D	Specimen			
41	Specimen35	49 AMP.M	007SP040.D	Specimen			
42				Keyword		NewBatch	
43	Blank02	50 AMP.M	007BL041.D	Blank			
44	CutOff02	51 AMP.M	007CU042.D	Cal	CU		
45	Negative02	52 AMP.M	007NE043.D	Negative			
46	Low Control02	53 AMP.M	007CO044.D	Control	LOW		
47	Med Control02	54 AMP.M	007CO045.D	Control	MED		

Step 3: Specify reports

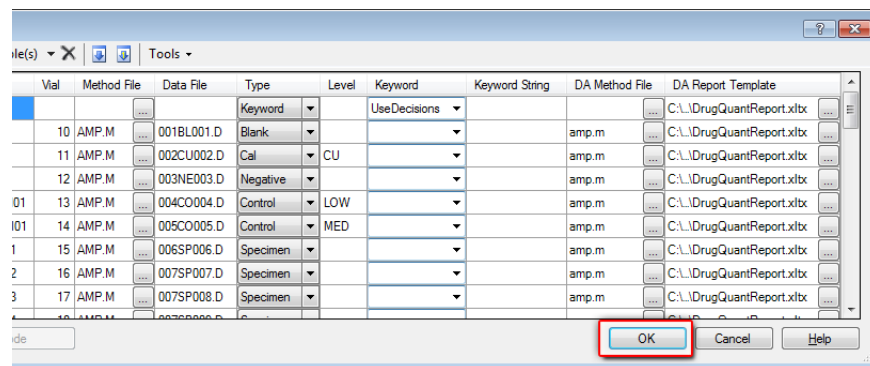
- In the **DA Report Template** column, click the browse button and select a report template.

This template is used to generate a report automatically with each sample.



Step 4: Save the Sequence Table

- Once you have completed entering all your samples in the Sequence Table, click **OK** to close the dialog.



- Save the completed Sequence Table.



Step 5: Run the Sequence

1. Load a previously created sequence table containing your batches of specimens.
2. Click the view icon to view the currently loaded sequence table and edit if needed.
3. With your sequence loaded, click the Start Sequence icon to display the Start Sequence dialog.
4. Notice that the name of the current sequence is displayed in the title bar.

For this Workflow, we are acquiring data, so we selected the **Full Method** option in the **Method Sections to Run** group box.

5. For the **Data File Directory** enter the path where you would like the batch directories to be created. Each automatically created batch directory will contain the acquired data files and a copy of the method used to acquire the data.
6. Click **Run Sequence** to begin processing all samples in this sequence.

In the previous sections you learned how to create a sequence table containing multiple batches of specimens. Once your sequence table has been set up, and your vials are loaded into the sample tray, you may begin to process your samples as described here.



Start Sequence DrugSpecimen.sequence.xml Last Modified: Wed May 22 11:03:11 2013

Method Sections to Run

Full Method

Reprocessing Only

Sequence Barcode Options

Disable barcode for this sequence.

On mismatch, inject anyway.

On mismatch, don't inject, continue the sequence.

On mismatch, don't inject, stop the sequence.

Overwrite Existing Data Files

Sequence Comment: _____

Operator Name: TWI\wtj

Data File Directory: C:\Quant\Batches\demo6\ Browse...

Pre-Sequence Macros/Commands

Acquisition: _____ Browse...

Data Analysis: _____ Browse...

Post-Sequence Macros/Commands

Acquisition: _____ Browse...

Data Analysis: _____ Browse...

Run the method as specified in the Method Information dialog box

Run Sequence OK Cancel Help

As each sample is analyzed, the results are compared with the parameters set in the method. If the results are outside the specified criteria, MassHunter GCMS Acquisition will continue the process based on the parameters specified in the Intelligent Sequencing portion of your method.

6

Update the Calibration

Introduction 76

Step 1: Create the Sequence Table 76

Step 2: Prepare the calibration samples 80

Step 3: Prepare the cutoff calibration sample 80

Step 4: Load the calibration sample vials in the ALS 80

Step 5: Run the Calibration Sequence 81

Update the Calibration

Create a
Sequence

Prepare
Samples

Run Samples



Agilent Technologies

Introduction

Step 1: Create the Sequence Table

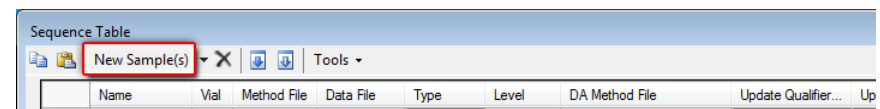
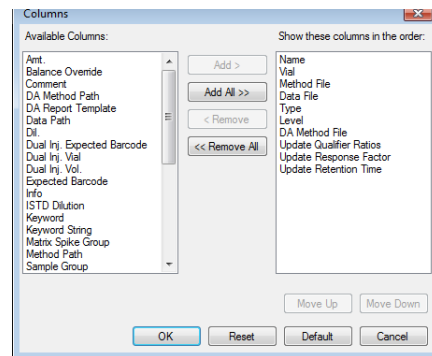
1. In the Instrument Control window, click the Edit Sequence icon to open the STE.
2. From the **Tools** menu select **Add/Remove Columns** and edit your columns to be similar to those shown here, then click **OK**.
3. Click the **New sample(s)** menu three times to add three blank lines to the sequence table.
4. The example shown here illustrates a completed table. We will describe how it was created in the following steps.

To begin, in the **Name** column, enter a sample name for each sample.

As required by governing regulations, the calibration curve stored in a method must be updated when a specified time has elapsed or when a quality control sample indicates an unacceptable deviation from the stored calibration curve.

This chapter describes how the Sequence Table Editor (STE) in MassHunter GCMS Acquisition can be used to automate this process.

In the following steps we will describe how you can modify this table to define your calibration samples.



	Name	Vial	Method File	Data File	Type	Level	DA Method File	Update Qualifier...
1	Level_100	11	AMP.M	AmpCal001	Cal	L100	Amp.quantmethod.xml	Replace
2	Cutoff Calibrator	12	AMP.M	CU002.D	Cal	CU	Amp.quantmethod.xml	Replace
3	Level_500	13	AMP.M	AmpCal002	Cal	L500	Amp.quantmethod.xml	Replace
4	Level_900	14	AMP.M	AmpCal003	Cal	L900	Amp.quantmethod.xml	Replace
5	Control_LOW	15	AMP.M	CTRL_LOW	Control	QC_LOW	Amp.quantmethod.xml	Replace
6	Control_MED	16	AMP.M	CTRL_MED	Control	QC_MED	Amp.quantmethod.xml	Replace

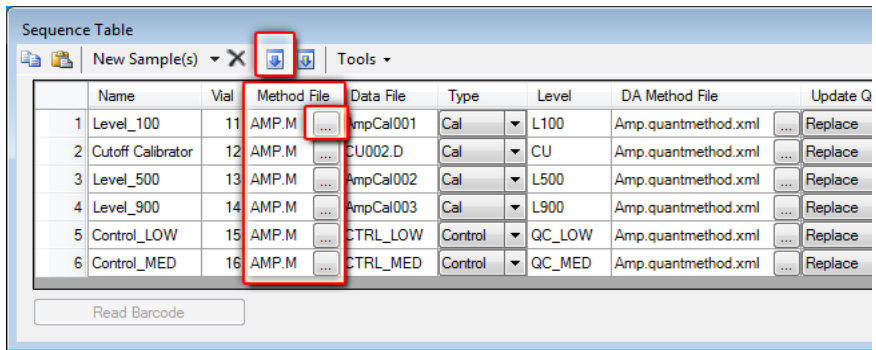
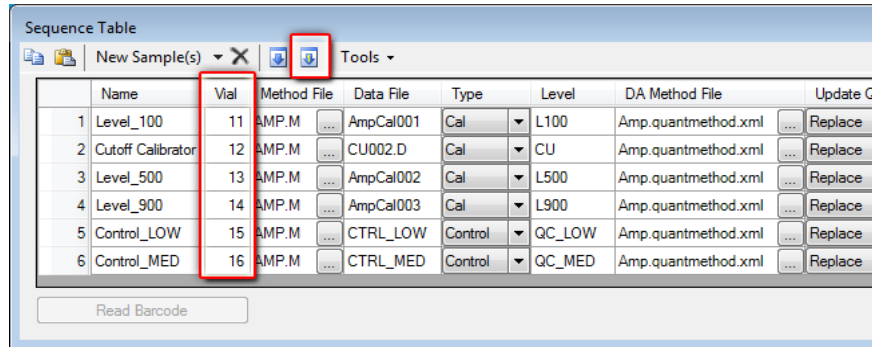
- Lines 1 through 4 are the 4 concentrations (**Levels**) required for a linear 4 point calibration curve for both Amphetamine and Methamphetamine. Line 2 is also a cutoff calibrator for these two drugs of abuse.
- Lines 5 and 6 are the control samples in our example.

5. In the **Vial** column, enter the location for the first sample (11 here), then click the increment icon (highlighted here), to automatically fill-down the vial numbers, (12 through 16 shown here).

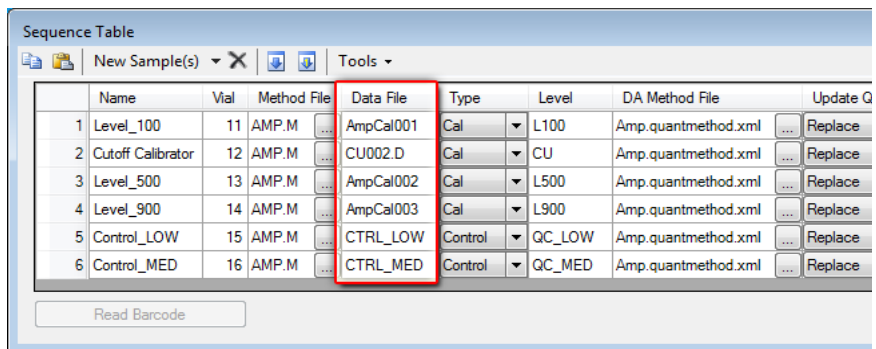
6. Click the browse icon next to the **Method File** name and select the method to use for each sample.

With your cursor on the first line, click the copy icon to copy this method name to each line in the table.

7. In the **Data File** column, enter a destination file name for the data acquired in the acquisition run.

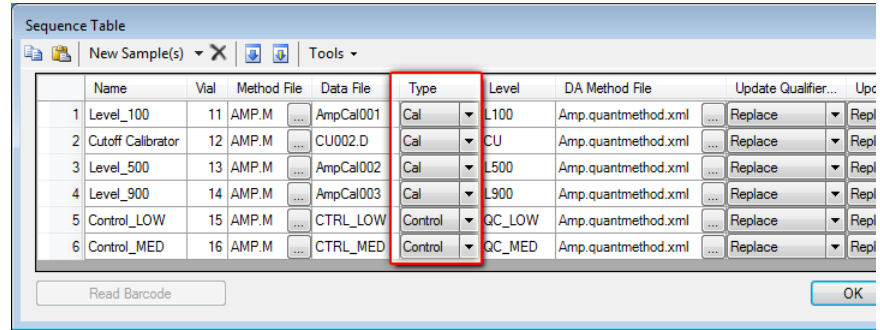


In this case, the instrument method parameters are stored in the AMP.M method, so that is the method to be used for each sample.



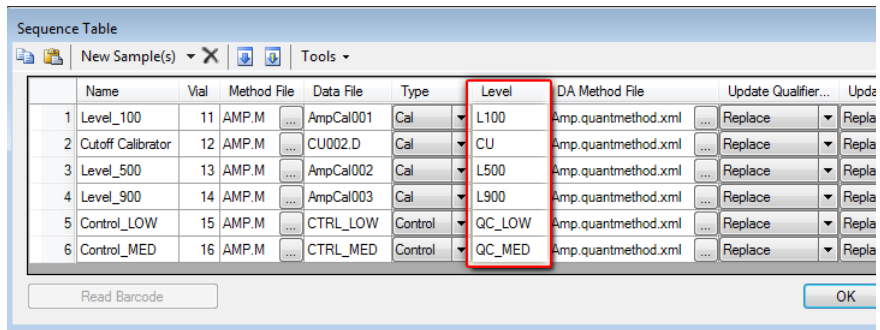
In this case the data will go into data files named AmpCal001, CU002, AmpCal002, AmpCal003, CTRL_LOW, and CTRL_MED.

8. In the **Type** column, select **CAL** from the dropdown list for lines 1 through 4 and select **Control** for lines 5 and 6.

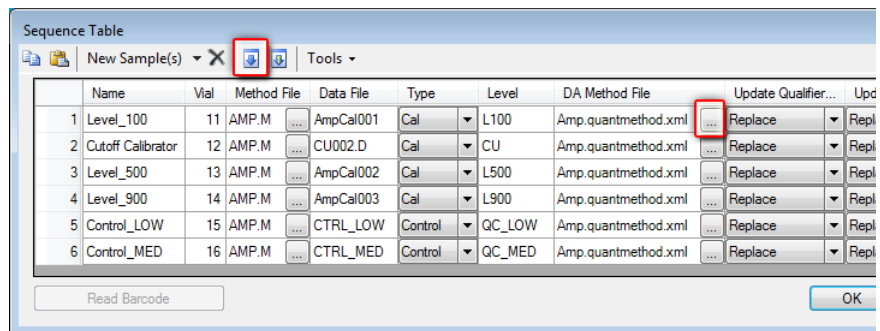


9. In the **Level** column, enter a level ID for each **CAL** and **Control** type sample.

This label must use *exactly the same name* as the corresponding concentration level ID in MassHunter Quantitative Analysis method. Capitalization is (is not) important, i.e. CU is not Cu.



10. In the **DA Method File** column, click the browse button on line 1, and select the MassHunter Quantitative Analysis method created in the last chapter. With the cursor still in this cell, click the copy icon to copy this method to all other samples in the table.



11. In the **Update Qualifier**, **Update Response**, and **Update Retention** columns, on line 1 select **Replace** from each dropdown. Copy this parameter to lines 2 through 6 for these three columns.

Vial	Method File	Data File	Type	Level	DA Method File	Update Qualifier...	Update Respon...	Update Retentio...
11	AMP.M	AmpCal001	Cal	L100	Amp.quantmethod.xml	Replace	Replace	Replace
12	AMP.M	CU002.D	Cal	CU	Amp.quantmethod.xml	Replace	Replace	Replace
13	AMP.M	AmpCal002	Cal	L500	Amp.quantmethod.xml	Replace	Replace	Replace
14	AMP.M	AmpCal003	Cal	L900	Amp.quantmethod.xml	Replace	Replace	Replace
15	AMP.M	CTRL_LOW	Control	QC_LOW	Amp.quantmethod.xml	Replace	Replace	Replace
16	AMP.M	CTRL_MED	Control	QC_MED	Amp.quantmethod.xml	Replace	Replace	Replace

The drop-down selections in these columns allow you to:

- **Replace** the existing parameter in the Quantitative Analysis method with the results from that sample analysis
- **Average** the result with the parameter in the Quantitative Analysis method
- **No Update** to analyze the CAL sample without updating the parameters in the Quantitative Analysis method

12. Click **OK** to close the sequence table.

Name	Vial	Method File	Data File	Type	Level	DA Method File	Update Qualifier...	Update Retention...
1 Level_100	11	AMP.M	AmpCal001	Cal	L100	Amp.quantmethod.xml	Replace	Replace
2 Cutoff Calibrator	12	AMP.M	CU002.D	Cal	CU	Amp.quantmethod.xml	Replace	Replace
3 Level_500	13	AMP.M	AmpCal002	Cal	L500	Amp.quantmethod.xml	Replace	Replace
4 Level_900	14	AMP.M	AmpCal003	Cal	L900	Amp.quantmethod.xml	Replace	Replace
5 Control_LOW	15	AMP.M	CTRL_LOW	Control	QC_LOW	Amp.quantmethod.xml	Replace	Replace
6 Control_MED	16	AMP.M	CTRL_MED	Control	QC_MED	Amp.quantmethod.xml	Replace	Replace

13. Save this sequence as a new name. Enter a new name for the sequence and save it to your instrument directory sequence folder.



Step 2: Prepare the calibration samples

Prepare calibration samples for creating a calibration curve of each drug of abuse compound in the MassHunter Quantitative Analysis method.

Each calibration sample should contain the drug of abuse compounds in concentrations that cover the expected linear range of each drug's calibration curve.

Each sample of a specified concentration should be spiked with those ISTD compounds referenced by each drug of abuse contained in the sample. In our example we are using 250 ng/mL of each ISTD in each of our CAL standards.

Our example uses a 4 point calibration requiring 4 concentrations of the compounds in individual vials. Here those concentrations are 100, 400, and 900 ng/mL for both Amphetamine and MethAmphetamine. The cutoff calibration sample contains the drugs and ISTD's for the 250 ng/mL level and is the second point on the calibration curve.

Step 3: Prepare the cutoff calibration sample

Prepare a cutoff calibration sample containing each drug of abuse in the concentration level required by the governing regulation. In our example the cutoff sample contains 250 ng/mL of both Amphetamine and Methamphetamine.

Spike the sample with those ISTD compounds as done with the other calibration samples.

Step 4: Load the calibration sample vials in the ALS

Once you have created your Sequence Table, place the vials into the sample tray and run the batch.

This example assumes use of an auto liquid sampler (ALS) which would require the 4 sample vials be placed in the sampler tray vial locations specified in the STE.

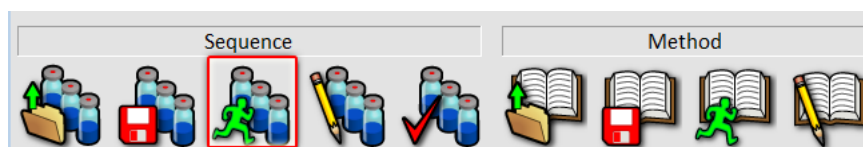
Step 5: Run the Calibration Sequence

1. With the newly created sequence loaded, click the **Run Sequence** icon.

2. On the Start Sequence dialog, select **Full Method** and enter a batch subdirectory for your acquired data files.

When the run is completed, the data files will be stored in the **Data File Directory**.

3. Click **Run Sequence** to begin the automated acquisition.
4. When the run is complete, in MassHunter Quantitative analysis **open this Batch** of 6 samples with the updated method and review the Calibration curve for linearity.
5. Enter the Quant Method Editor and examine these updated results in the method table.
6. Exit MassHunter Quantitative Analysis.



Start Sequence AmpCalibration.sequence.xml Last Modified: Wed May 15 15:59:43 2013

Method Sections to Run

Full Method

Reprocessing Only

Sequence Barcode Options

Disable barcode for this sequence.

On mismatch, inject anyway.

On mismatch, don't inject; continue the sequence.

On mismatch, don't inject; stop the sequence.

Overwrite Existing Data Files

Sequence Comment: _____

Operator Name: TWI\wit

Data File Directory: C:\Quant\Batches\demo2\

Pre-Sequence Macros/Commands

Acquisition: _____

Data Analysis: _____

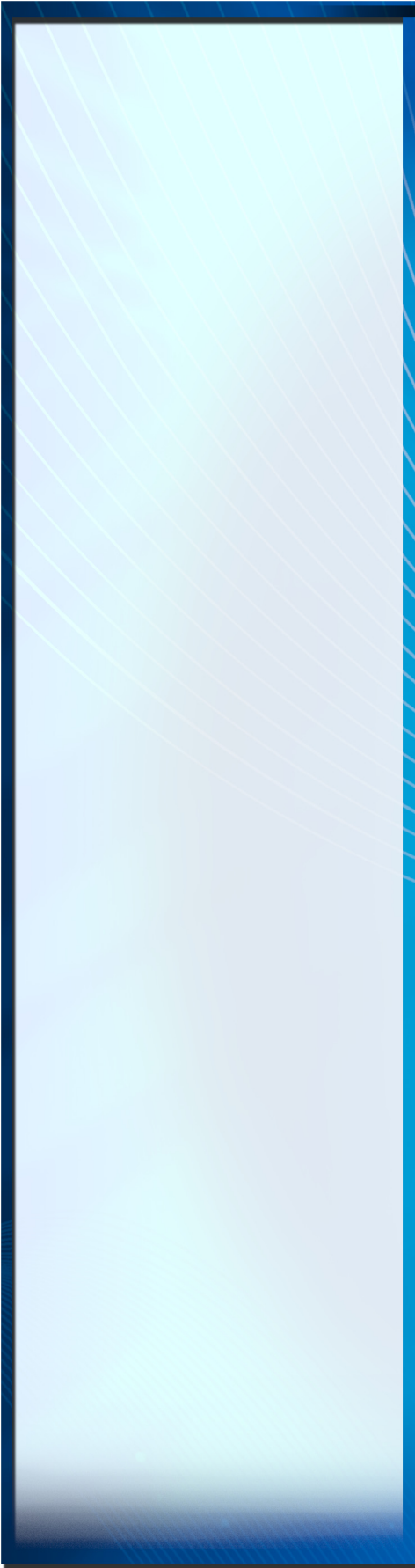
Post-Sequence Macros/Commands

Acquisition: _____

Data Analysis: _____

Run the method as specified in the Method Information dialog box

The data for a sample is acquired and analyzed before the next sample in the table is processed. The quantitative method is updated by replacing the qualifier ratios, detector response, and retention time of the quantitative ion with the actual results from the analysis of the sample. This is done for the level specified for this sample in the Sequence table.





© Agilent Technologies, Inc.

Printed in USA, August 2013