

Axiom™ Analysis Suite 3.1

USER GUIDE

Publication Number 703307

Revision 5

The information in this guide is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Legal entity

Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

TRADEMARKS

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2017 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■	CHAPTER 1 Introduction	9
	Overview	9
	Software and Hardware Requirements	9
	Sample Data Size Estimates and Required Disk Space	10
	Installation Instructions	11
	Starting Axiom Analysis Suite	12
	Using the Preferences Window Tab	14
	Changing the Default Library Folder/Path	15
	Setting Up Proxy Server Access	15
	Updating NetAffx Library/Annotations	16
	Enabling/Disabling Check for Library File Updates at Start Up	17
	Installing Custom Array Library Files	17
	Uninstalling	18
	Windows 7	18
	Windows 10	18
■	CHAPTER 2 Performing an Analysis	19
	Setting Up an Analysis	19
	Selecting a Mode (Workflow)	19
	Selecting an Array Type	20
	Importing CEL Files	20
	Importing CEL Files by Text	21
	Removing Selected CEL Files	21
	Setting Up an Analysis Configuration	22
	Selecting an Analysis Configuration	22
	Using the Analysis Settings Fields	24
	Sample QC Fields	24
	Genotyping Fields	24
	Saving your Analysis Configuration	26
	Modifying an Existing Analysis Configuration	27
	Setting Up Threshold Settings	28
	Customizing Thresholds	28
	Sample QC	29
	SNP QC	29
	Assigning an Output Folder Path	31
	Assigning a New Output Folder Path	31
	Adding Sub-Folders	31
	Assigning a Batch Name	31

Running your Analysis	32
Using the Dashboard Window Tab	34
Open Selected Result(s)	35
Remove Selected Result(s)	35
Viewing the Results Folder in Windows Explorer	35
Browsing For Existing Analysis Results	36
Browsing for Existing Suitcases	37
Opening an Analysis from Windows Explorer	37

■ **CHAPTER 3 The Viewer: Summary Window and Sample Table . . 38**

Viewing Options	39
Split-Screen Options	39
Changing a Tab Window to a Full Screen Windows	41
Adjusting the Window Size	42
Summary Window/Tab	43
Data Analysis Summary	43
Viewing the Plate Barcode Table Details	44
Sample Table	45
Importing Sample Attributes	46
Column Headers	46
Rearranging Columns	47
Sorting Columns	47
Single-Click Sorting Method	47
Hiding the Column	47
Filtering Column Data	48
Adding Filters (Method 1)	48
Text-based Columns	48
Numeric Data Columns	49
Showing Filtered Data Only	50
Clearing an Individual Filter	51
Clearing All Current Filters	52
Adding Filters (Method 2)	52
Copying Column Data	55
Setting User Colors	55
Assigning a Color to a Sample	55
Importing Assigned Colors	56
Viewing User Colors in the Cluster Graph	58
Removing an Assigned User Color	59
Searching Keywords	60
Box Plots	61
Viewing the Default Box Plots	61
Changing the Box Plot's Scale Setting Ranges	62
Adding a New Box Plot	62
Reading Box Plot Percentiles	63
Saving the Current Box Plot View	63

Scatter Plot	64
Viewing the Default Scatter Plot	64
Changing the Scatter Plot's Setting Ranges	65
Adding a New Scatter Plot and Selecting its X and Y Properties	65
Customizing Color By Settings	67
Saving the Current Scatter Plot View	68
Plate Views	69
Viewing the Default Plate Views	69
Adding a New Plate View Metric	70
Customizing your Plate View Settings	71
Saving the Current Plate View	71
Concordance Checks	72
Running a Concordance Check	72
Comparing All Combinations	72
Comparing to Reference	74
Reanalyzing Samples	74
■ CHAPTER 4 The Viewer: SNP Summary Table and Cluster Plot .	76
SNP Summary Table	76
Using the SNP Summary Table	78
Setting your SNP Summary Table View	78
Adding and Removing Table Columns	78
Selecting Annotations	79
Saving your Table Column View	80
Copying Selected Row(s)	80
Copying Selected Cell(s)	80
Changing or Reverting Genotype Calls	81
Changing Genotype Calls by Text File	81
Reverting Selected SNPs for all Samples	81
Reverting all SNPs for all Samples	81
Reanalyzing your SNP Summary Table Data	82
Regenerate SNP Metrics	82
Running OTV Caller	85
Managing your SNP List	86
Saving your Current SNP List	86
Exporting your SNP List	86
Importing a SNP List	87
Using your Saved SNP List	87
Displaying SNPs in a SNP list	87
Displaying SNPs that are not in your SNP List	88

Cluster Plot	89
Using the Cluster Plot	91
Displaying a SNP Cluster Plot that Corresponds with a SNP	91
Setting New Scale Setting Ranges	92
Customizing Color By Settings	92
Selecting Multiple Samples in a Cluster Plot	94
Changing a Sample's Call for a Single SNP	95
Reverting a Single Call	95
Reverting Multiple Calls	95
Displaying Cluster Model Data	96
Saving the Current Cluster Plot View	96
■ CHAPTER 5 Allele Translation	97
About Translations	97
Performing Allele Translation	97
Allele Translation Options	99
Translation Reports	101
Comprehensive and Summary Translation Report	102
Summary Translation Report	102
Phenotype Translation Report	103
Phenotype Report	103
Opening Translation Report in MS Excel	104
Available Report Fields and Descriptions	105
Array Tracking	105
Gene-specific	105
Marker-specific	108
Tracking Edited Genotype Calls	109
Uncalled Report	110
■ CHAPTER 6 Exporting	111
Using the Sample Table Export Options	111
Using the SNP Summary Table Export Options	111
Exporting the Current Table	112
Exporting All Data	112
Exporting Signature SNPs	112
Exporting Genotyping Data	112
Result Output Formats	113
Call Output Formats	114
Exported Data Selections	114
Input and Output Files	114

SNP List Filter (Optional)	115
Output Location (Required)	115
Output Name (Required)	115
Changing the SNP Identifier	115
Changing the Current Annotation File (Optional)	116
Adding and Removing Annotation Columns	116
Exporting Cluster Plots to PDF	117
■ CHAPTER 7 External Tools	119
Axiom CNV Tool 1.1	120
Council on Dairy Cattle Breeding (CDCB) Export Tool	120
Axiom Long Format Export Tool	121
■ APPENDIX A Predefined Region CN Analysis Variations	122
Predefined Copy Number Analyses	122
With Predefined CN Genotyping	122
With CN-aware Genotyping	123
Genotyping with CN Analysis	124
Setting Up a Genotyping with CN Analysis	124
Selecting a Mode (Workflow)	124
Importing CEL Files	124
Analysis Settings with Copy Number Options	127
Sample QC	127
Genotyping	127
Threshold Configurations with CN Options	129
Sample QC	129
CN QC	129
SNP QC	129
Assigning an Output Folder Path	129
Assigning a Batch Name	129
Running your Genotyping with CN Analysis	130
Viewing your Genotyping and CN Results	130
Summary Report	131
Sample Table	133
SNP Summary Table	134
CN Summary Table and CN Region Plot	135
CN Summary Table (Overview)	135
CN Region Plot (Overview)	136
Overview and Use of the Best Practices Workflow	137

■ APPENDIX B About Allele Translation.....	139
Overview	139
Gene Table Layout for Haplotyping	139
Biological Annotations in Translation Reports	141
Impact of Phase Ambiguity in Haplotyping	142
Diplotype to Phenotype Translation	144
Creating a Custom Metabolizer Library File	145
Metabolizer Library File Format	146
Reference Databases Used in Translation Data Curation	148
■ APPENDIX C Definitions	149
Threshold Names	149
Sample Table	154
Concordance Columns	158
Annotations and Columns	159
SNP Summary Table Definitions	162



Introduction

Overview

Axiom Analysis Suite (AxAS) enables you to perform the following functions:

- Run QC and Genotyping Algorithms.
- View QC Data within tables and graphs at a Sample and/or SNP level.
- View Cluster Graphs with the ability to change calls and/or highlight by attribute.
- Export your Data.

Software and Hardware Requirements

64-bit Operating System	Speed	Memory (RAM)	Available Disk Space ¹	Web Browser
Microsoft Windows® 7 (64 bit) Professional with Service Pack 1	2.83 GHz Intel Pentium Quad Core Processor	16 GB	150 GB HD + data storage Refer to table below.	Internet Explorer 8.0 and above
Microsoft Windows 10 (64 bit) Professional	2.83 GHz Intel Pentium Quad Core Processor	16 GB	150 GB HD + data storage Refer to table below.	Internet Explorer 8.0 and above

¹Minimum storage requirements are for a single run. Total storage space should include additional space for data storage of input and output files from current and previously completed analyses. In addition, you must have a minimum of 5GB of free space on your C: drive to run an analysis.

Sample Data Size Estimates and Required Disk Space

Before using AxAS, make sure you have enough disk space. See the table below for size estimates. Note: The estimates shown include the contents of the batch name folder.²

# of Markers	Storage Type ³	50 samples	100 samples	500 samples	1000 samples	5000 samples
50K	Input	1.33 GB	2.66 GB	13.3 GB	26.6 GB	133 GB
	Output	158 MB	286 MB	1.27 GB	2.51 GB	12.4 GB
	Total	1.49 GB	2.95 GB	14.57 GB	29.11 GB	145.4 GB
500K	Input	1.33 GB	2.66 GB	13.3 GB	26.6 GB	133 GB
	Output	1.53 GB	2.77 GB	12.6 GB	25.0 GB	124 GB
	Total	2.86 GB	5.43 GB	25.9 GB	51.6 GB	257 GB
850K	Input	1.33 GB	2.66 GB	13.3 GB	26.6 GB	133 GB
	Output	2.59 GB	4.69 GB	21.4 GB	42.4 GB	209 GB
	Total	3.92 GB	7.35 GB	34.7 GB	69.0 GB	342 GB

²A batch name folder is auto-generated during the analysis process. This folder includes all the necessary files needed to view your analysis results in the Viewer.

³Input is the storage size required for CEL files to be analyzed. Output is the storage size required for analysis results files.

Installation Instructions

1. Go to thermofisher.com, then navigate to the following location:
Applications and Techniques > Life Sciences > Microarray Analysis > Software > Axiom Analysis Suite

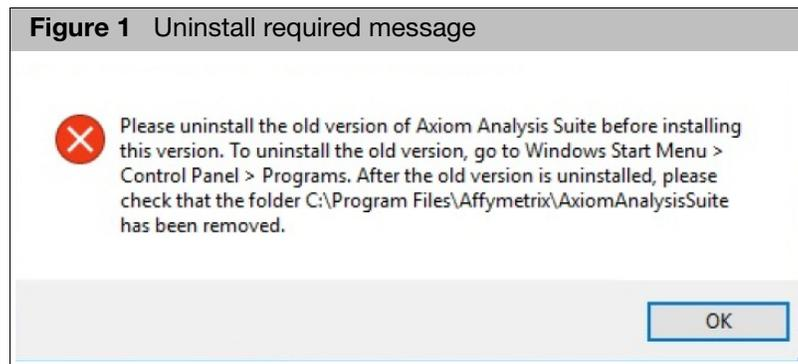
Or

Click on this link:

<https://www.thermofisher.com/us/en/home/life-science/microarray-analysis/microarray-analysis-instruments-software-services/microarray-analysis-software/axiom-analysis-suite.html>

2. Locate and download the zipped Axiom Analysis Suite software package.
3. Unzip the file, then double-click **AxiomAnalysisSuiteSetup.exe**.
4. Follow the on-screen instructions to complete the installation.

If your system has a previous version installed, the following message appears:
(Figure 1)

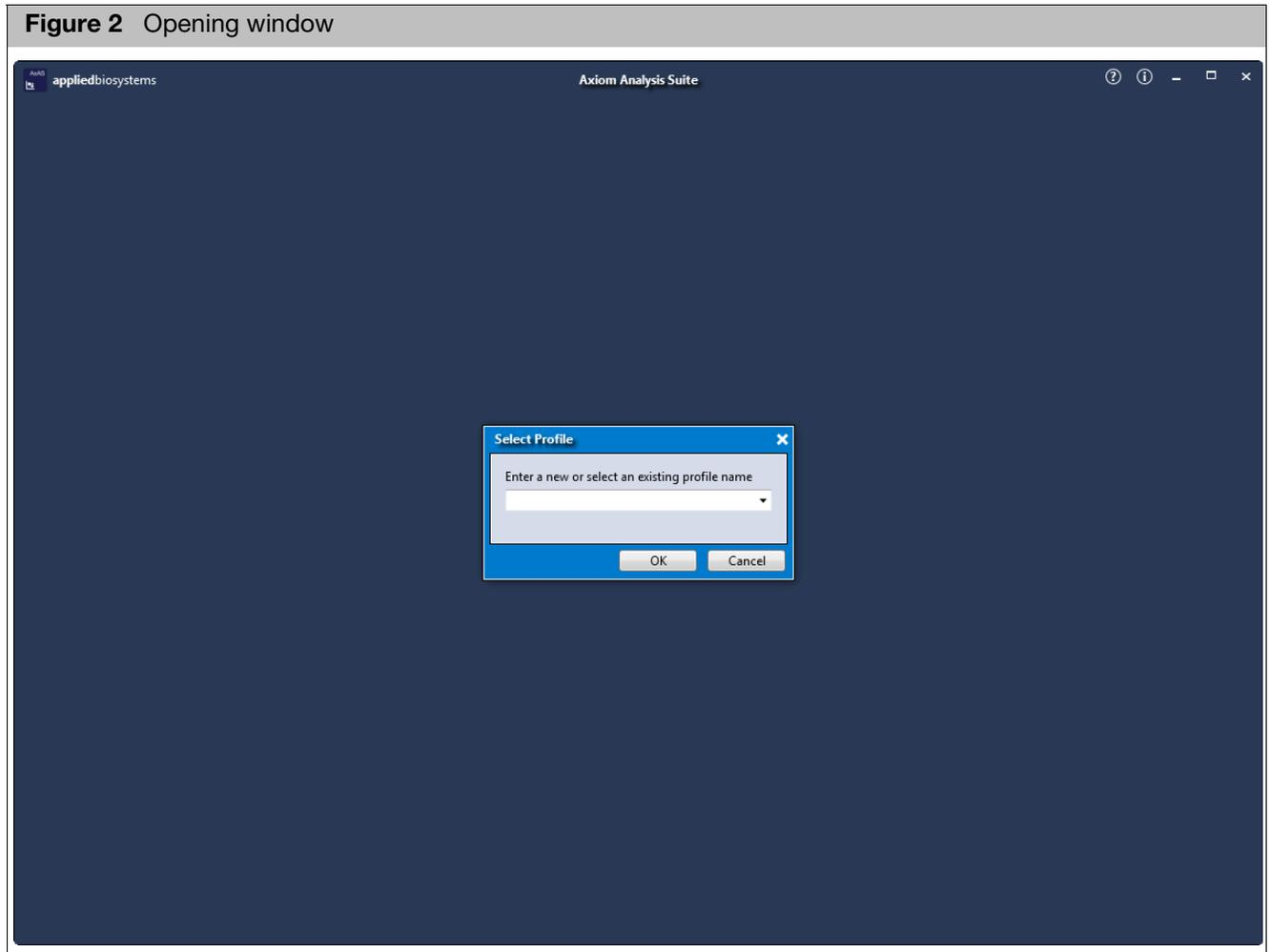


Acknowledge the message, click **OK**, then go to "Uninstalling" on page 18.

Starting Axiom Analysis Suite

1. Double-click on the Axiom Analysis Suite Desktop shortcut or click **Start > All Programs > Thermo Fisher Scientific > Axiom Analysis Suite**.

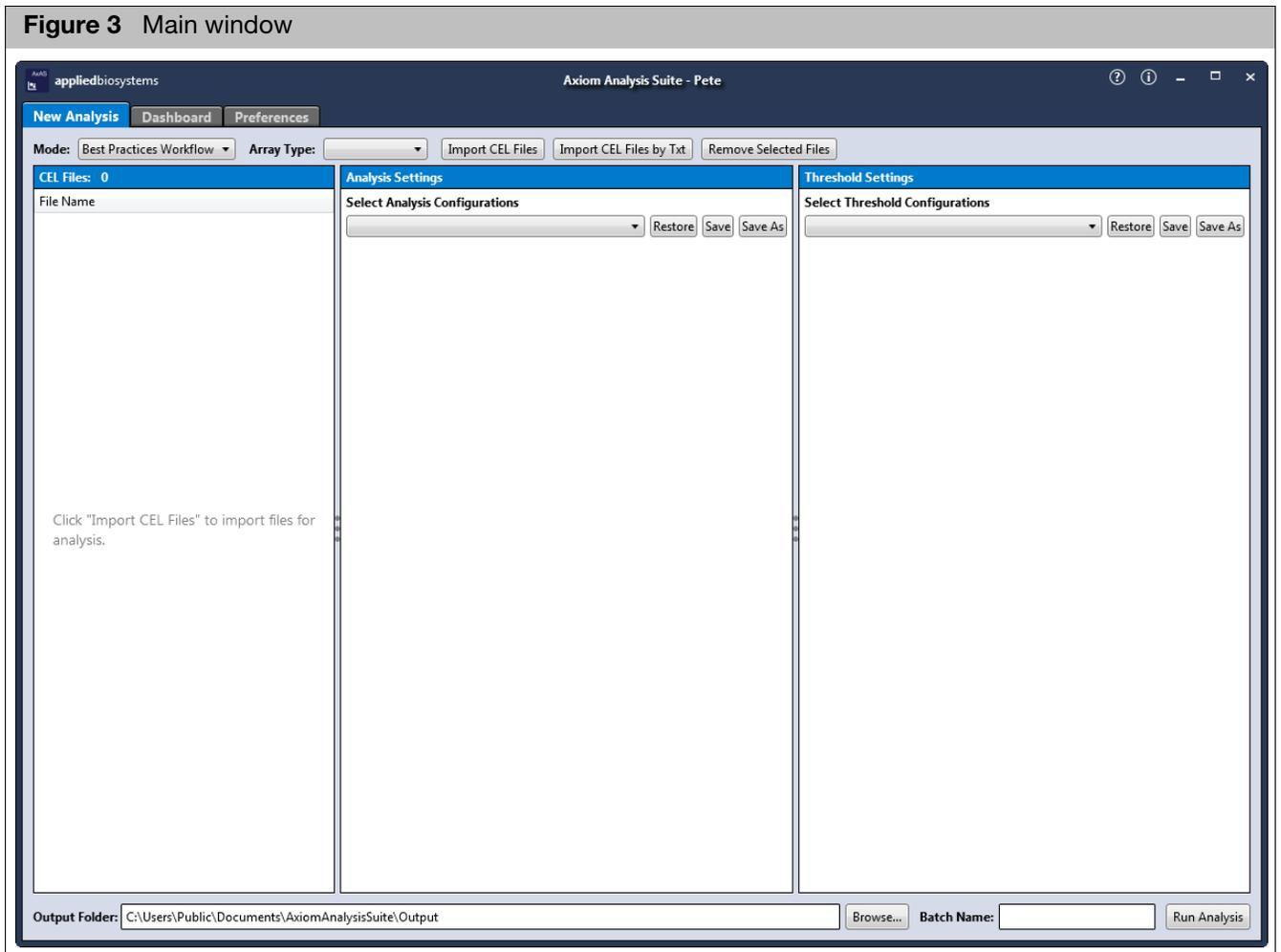
The following window appears: (Figure 2)



2. Enter a new profile name or click the down-arrow to select an existing profile name.
3. Click **OK**.

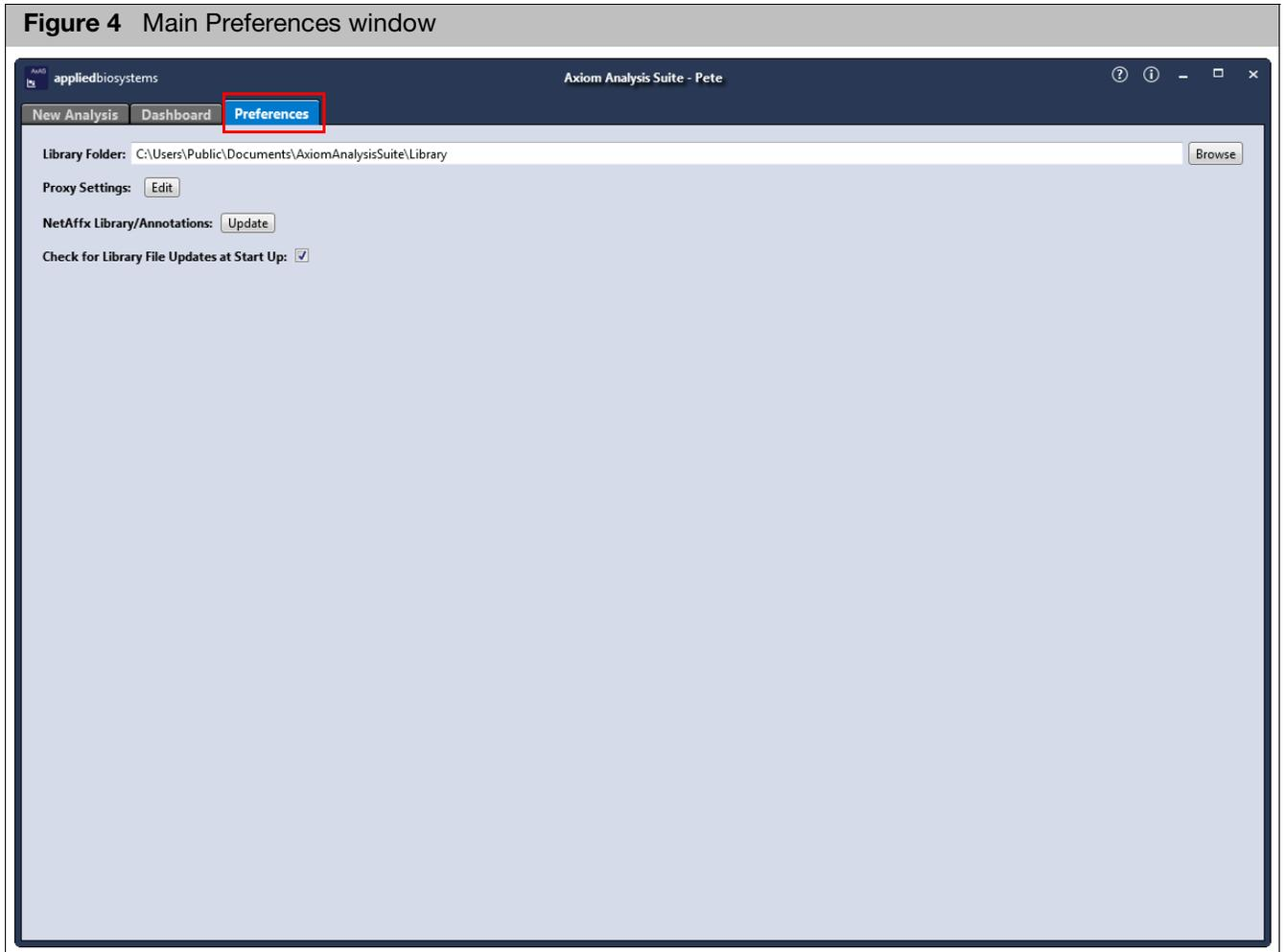
The following window appears: (Figure 3)

Figure 3 Main window



Using the Preferences Window Tab

Click the **Preferences** window tab (Figure 4) to setup or change a library path, edit Proxy settings, download or update Library/Annotation files.

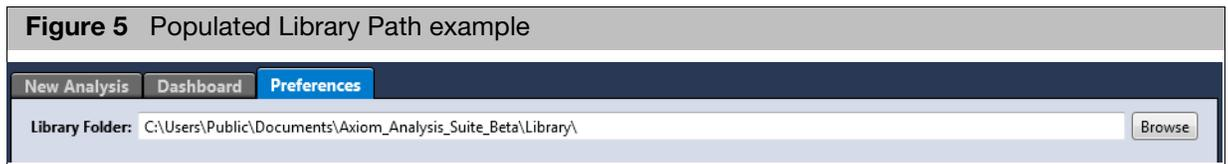


Changing the Default Library Folder/Path

✗ IMPORTANT! The library folder contains the library and annotation files required to run the Axiom Analysis Suite software.

Do the following to change the default Library folder/path:

1. Click **Browse** (right of library path field).
The **Select Library Folder** window appears.
2. Navigate to the new location you want the library folder to reside.
3. Click **New Folder**.
4. Rename the New Folder (as you normally would), then click **Select Folder**.
Your newly assigned Library folder is set and reflected in the Library Folder directory/path field, as shown in [Figure 5](#).

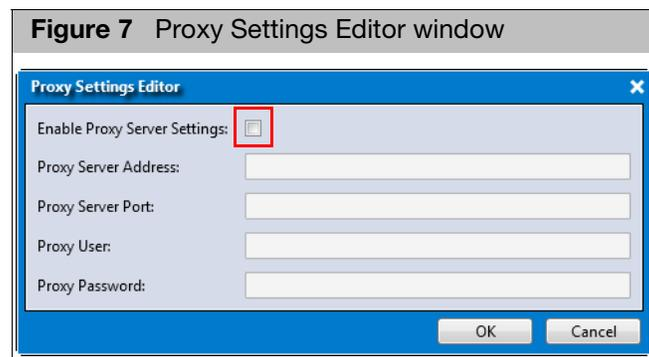


Setting Up Proxy Server Access

If your system has to pass through a Proxy Server before it can access the NetAffx server (Internet), click the **Edit** button. ([Figure 6](#))



The following window appears: ([Figure 7](#))



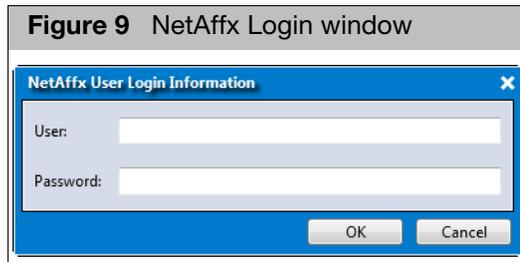
5. Click the **Enable Proxy Server Settings** check box ([Figure 7](#)), then contact your IT department for help with completing the required text fields.
6. Click **OK**.

Updating NetAffx Library/Annotations

1. Click on the **Update** button. (Figure 8)



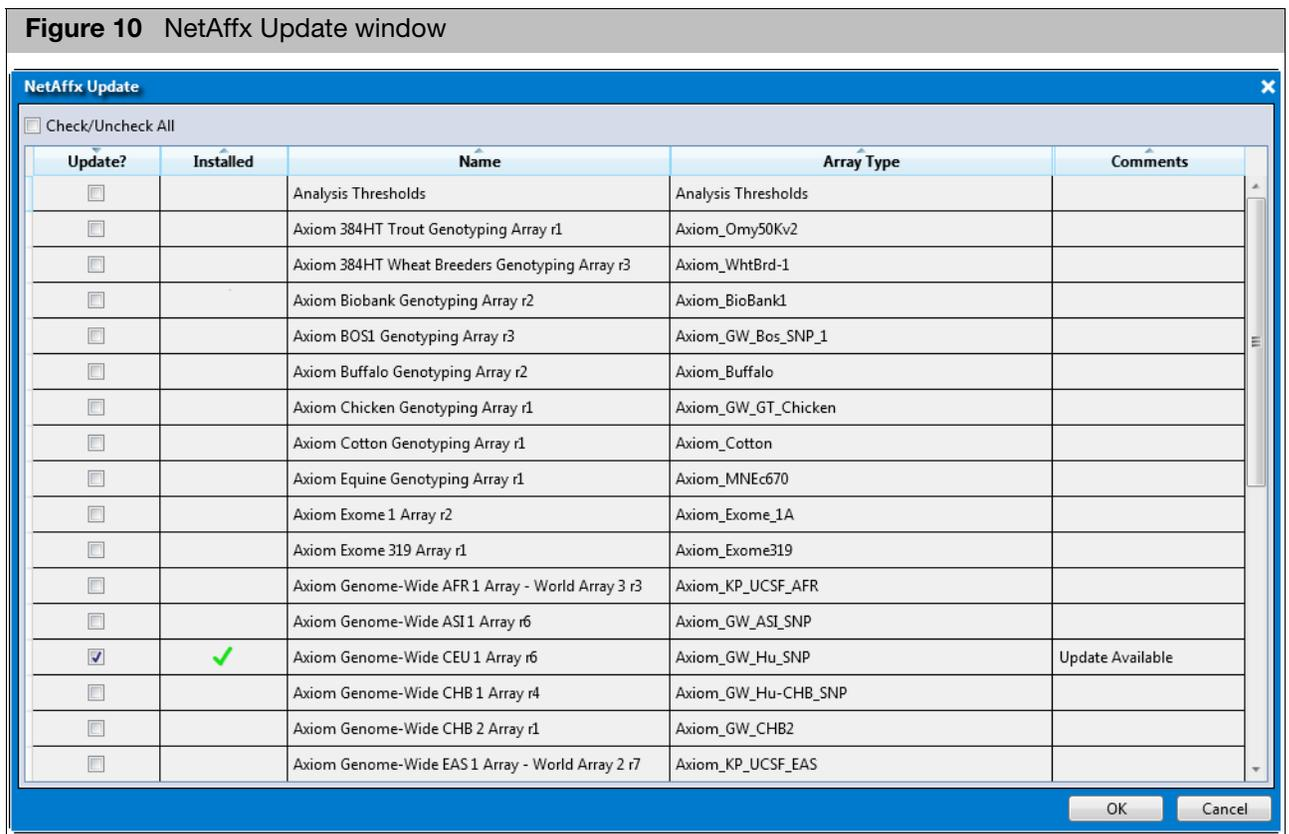
The following window appears: (Figure 9)



2. Enter your NetAffx account email and password, then click **OK** or go to www.netaffx.com and click **Register** to sign up.

Note: If you are unable to connect to NetAffx, make sure you have an active Internet connection, and/or correct Proxy Server settings.

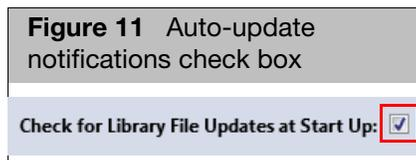
The NetAffx Update window appears. (Figure 10)



3. You must click the check box(es) that correspond with the type of CEL files you want to analyze.
Click the **Check/Uncheck All** check box to select/deselect all the listed check boxes.
4. Click **OK**.
An Installing Updates progress bar appears.

Enabling/Disabling Check for Library File Updates at Start Up

1. This check box (Figure 11) is checked by default to enable automatic Library File update alerts each time you launch the Axiom Analysis Suite application.
(Recommended)



Installing Custom Array Library Files

X IMPORTANT! Library files for custom designs must be manually installed.

1. Download the zip package provided to you by Affymetrix Bioinformatics Services.
2. Unzip the contents of the analysis library files into a single sub-folder within the library file folder.

For multi-species designs, each species should be in its own sub-folder. There should be no other folders within each sub-folder and all annotation information must be in the same location as the .CDF file.

Uninstalling

X IMPORTANT! Upgrade installations are NOT supported. You must uninstall the existing version of Axiom Analysis Suite BEFORE installing this new version. Administrative rights to the computer are required before you can uninstall the Axiom Analysis Suite software. For your convenience, no existing library files or user settings are removed during the uninstall process.

Windows 7

1. Click **Start > Control Panel**.
The Control Panel window appears.
2. Click the **View by** drop-down menu (upper-right), then click to select **Category**.
3. In the **Programs** category, click **Uninstall a program**.
The Programs and Features window appears.
4. Click to select **Axiom Analysis Suite**, then click **Uninstall**.
5. Follow the on-screen instructions.
6. After the uninstall process is complete, close the Programs and Features window.
7. Use Windows Explorer as you normally would to navigate to the directory:
C:\Program Files\Affymetrix
8. Verify that the **Axiom Analysis Suite** folder has been removed.
9. If the folder is present, double-click on it to open it.
10. Search for any files you want to keep, then move them to different (easily accessible) location.
11. Delete the **Axiom Analysis Suite** folder.
12. Close all open windows, then install the new version, as described in the "[Installation Instructions](#)" on page 11.

Windows 10

1. Click the Windows icon (bottom left corner).
2. Click **All apps > Windows System > Control Panel**.
The Control Panel window appears.
3. In the **Programs** category, click **Uninstall a program**.
The Programs and Features window appears.
4. Click to select **Axiom Analysis Suite**, then click **Uninstall**.
5. Follow the on-screen instructions.
6. After the uninstall process is complete, close all open windows.
7. Use Windows Explorer as you normally would to navigate to the directory:
C:\Program Files\Affymetrix
8. Verify that the **Axiom Analysis Suite** folder has been removed.
9. If the folder is present, double-click on it to open it.
10. Search for any files you want to keep, then move them to different (easily accessible) location.
11. Delete the **Axiom Analysis Suite** folder.
12. Close all open windows, then install the new version, as described in the "[Installation Instructions](#)" on page 11.

2

Performing an Analysis

After downloading the library and annotation files that match the array type of the CEL files you want to analyze, click the **New Analysis** tab.

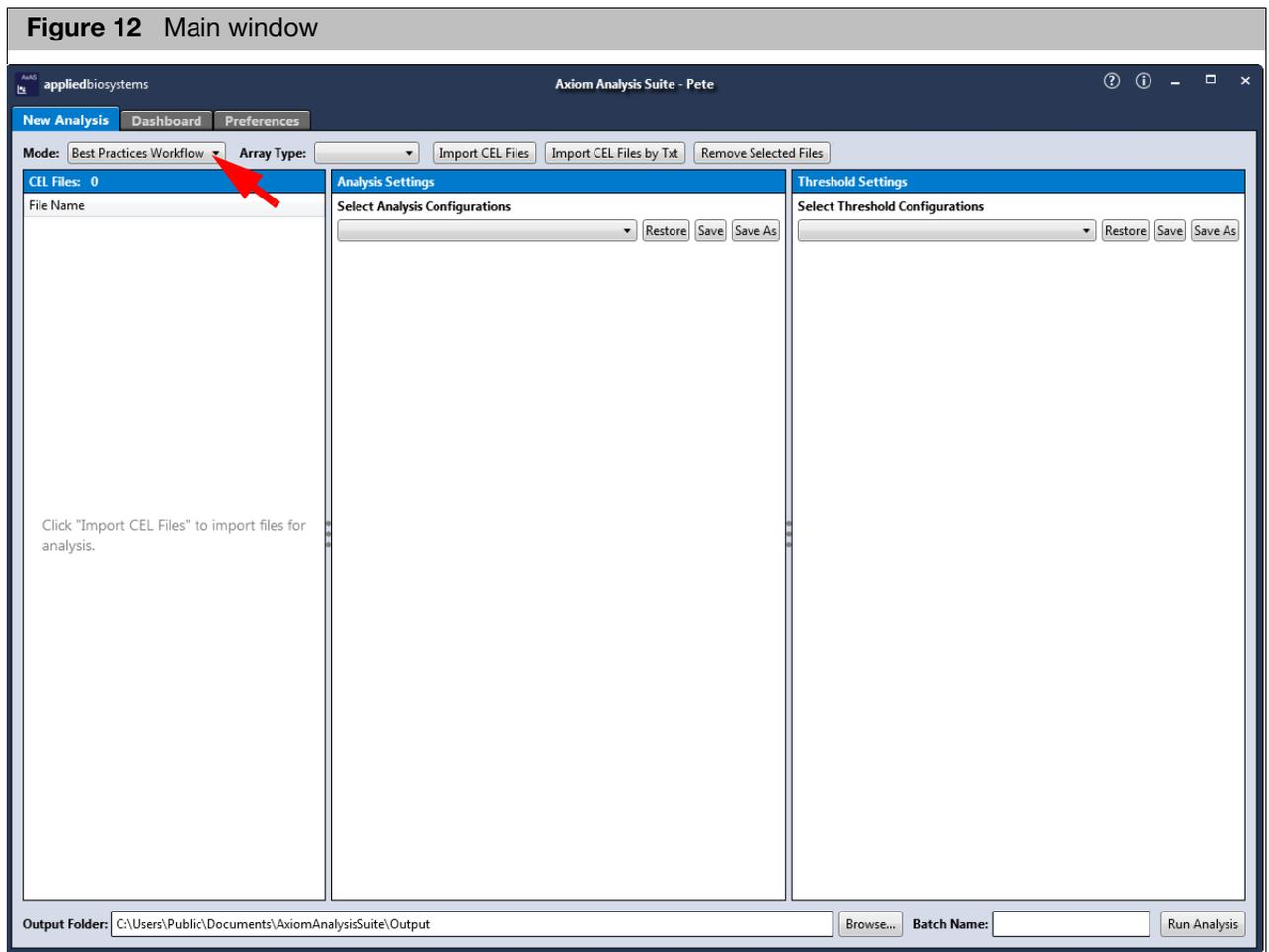
The **New Analysis** window and its three individual panes appear, as shown in [Figure 12](#).

Setting Up an Analysis

If you want to run and view CN-aware genotypes, go to [page 122](#). If not, continue to "Selecting a Mode (Workflow)".

Selecting a Mode (Workflow)

1. From the main Axiom Analysis Suite window tab, click the **Mode** drop-down. ([Figure 12](#))



2. Click to select the workflow you want to use.
 - **Best Practices Workflow (Default):** This workflow performs quality control analysis for samples and plates, genotypes those samples which pass the defined QC thresholds, and then categorizes the probe sets to identify those whose genotypes are recommended for statistical tests in downstream study. Details are available in the Axiom Genotyping Solution Data Analysis Guide (P/N 702961)
 - **Sample QC:** This workflow performs the quality control analysis for samples and plates. Note this workflow does not produce genotype calls for the passing samples.
 - **Genotyping:** This performs genotyping on the imported CEL files, regardless of the sample and plate QC metrics. Note: Including samples that do not pass defined QC thresholds may reduce the quality of the results for passing samples.
 - **Summary Only:** This workflow produces a summary of the intensities for the probe sets for use in copy number analysis tools. Note: Summary Only does not perform sample QC nor genotyping.

Selecting an Array Type

1. Click the **Array Type** drop-down to select the array type you want to use.

Importing CEL Files

1. Click **Import CEL Files**.
The **Add CEL Files** window appears.
2. Navigate to your CEL file location. Make sure the CEL Files you select coincide with the array type you selected earlier, otherwise a warning message appears.
3. Single-click on a CEL file or Ctrl click, Shift click, or press Ctrl A (to select multiple files).
4. Click **Open**.

The CEL Files pane populates and displays your selected CEL files. (Figure 13)

Figure 13 Populated CEL File pane example

CEL Files: 12
File Name
2877_AxiomGWHuSNP1_20101115_Blood_G01
2878_AxiomGWHuSNP1_20101115_Blood_H01
2879_AxiomGWHuSNP1_20101115_Blood_A02
2880_AxiomGWHuSNP1_20101115_Blood_B02
2881_AxiomGWHuSNP1_20101115_Blood_C02
2882_AxiomGWHuSNP1_20101115_Blood_D02

Importing CEL Files by Text

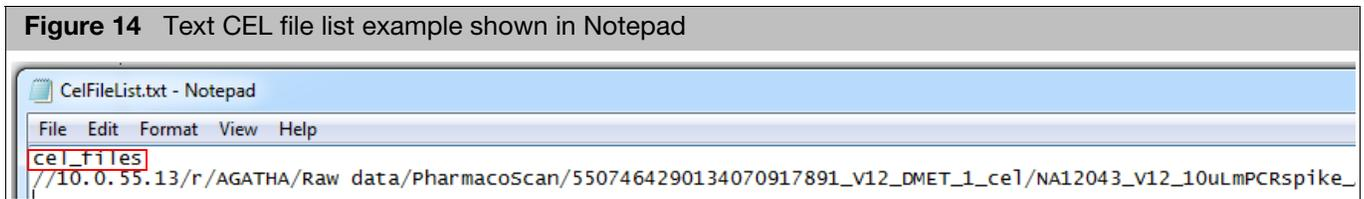
✗ IMPORTANT! The CEL files paths in a TXT file are not recognized (and will not import) if they begin with back slashes (e.g: `\\path\to\my\file.cel`).

Before importing, the two starting back slashes must be replaced with forward slashes (e.g: `//path\to\my\file.cel`).

Also, your `*.txt` list file must start with the header `cel_files`, as shown in [Figure 14](#).

1. Click **Import CEL Files by Txt**.
The **Import CEL Files by Txt** window appears.
2. Navigate to the `.txt` file that contains the list of CEL files you want to process.

Figure 14 Text CEL file list example shown in Notepad



Make sure the CEL Files you select coincide with the array type you selected earlier, otherwise a warning message appears.

3. Single-click on a CEL file or Ctrl click, Shift click, or press Ctrl A (to select multiple files).
4. Click **Open**.
Your CEL Files pane populates and displays each CEL file extracted from your selected text file.

Removing Selected CEL Files

Use this option to remove unwanted CEL files.

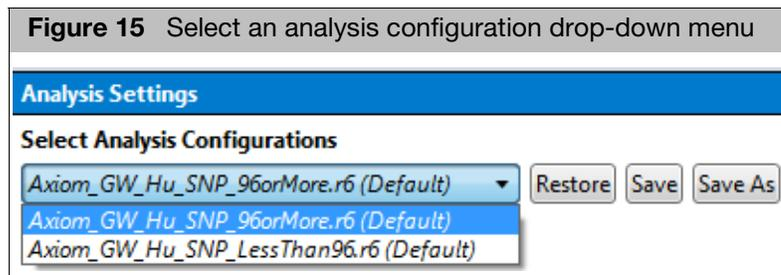
1. Single-click on a CEL file or Ctrl click, Shift click, or press Ctrl A (to select multiple files), then click **Remove Selected Files**.

Setting Up an Analysis Configuration

The Analysis Settings are populated based on the Mode (Workflow) chosen. For example, if Genotyping mode is selected, the Sample QC section of the Analysis Settings is hidden and only the Genotyping section is visible.

Selecting an Analysis Configuration

1. It is highly recommended you click the drop-down menu (Figure 15) and select the option that best matches the number of samples you want to analyze.
Note: The default configuration options displayed in the drop-down menu are based on your array type.



After selecting the appropriate default for the number of your samples, the Analysis Setting pane auto-populates, as shown in Figure 16.

Figure 16 Auto-populated Analysis Setting pane example

The screenshot displays the 'Analysis Settings' window with the following sections and fields:

- Select Analysis Configurations:** A dropdown menu showing 'Axiom_PMRA_96orMore.r1 (Default)', with 'Restore', 'Save', and 'Save As' buttons.
- Sample QC:**
 - Analysis File:** A dropdown menu showing 'Axiom_PMRA_96orMore_Step1.r1.appt-genotype-axiom.AxiomGT1.appt2'.
 - Prior Model File:** A text field containing 'Axiom_PMRA.r1.generic_prior' with browse and delete buttons.
 - SNP List File:** A text field containing 'Axiom_PMRA.r1.step1' with browse and delete buttons.
 - Gender File (optional):** An empty text field with browse and delete buttons.
 - Hints/Inbred File (optional):** Radio buttons for 'Inbred' and 'Hints' (selected), followed by an empty text field with browse and delete buttons.
- Genotyping:**
 - Analysis File:** A dropdown menu showing 'Axiom_PMRA.r1.appt-genotype-axiom.AxiomGT1.appt2'.
 - Prior Model File:** A text field containing 'Axiom_PMRA.r1.generic_prior' with browse and delete buttons.
 - Multi-allele Background Prior Model File:** A text field containing 'Axiom_PMRA.r1.generic_prior.mmb' with browse and delete buttons.
 - Multi-allele Pairwise Prior Model File:** A text field containing 'Axiom_PMRA.r1.generic_prior.mmp' with browse and delete buttons.
 - Multi-allele Prior Model File:** A text field containing 'Axiom_PMRA.r1.generic_prior.mm' with browse and delete buttons.
 - SNP List File:** A text field containing 'Axiom_PMRA.r1.step2' with browse and delete buttons.
 - Gender File (optional):** An empty text field with browse and delete buttons.
 - Hints/Inbred File (optional):** Radio buttons for 'Inbred' and 'Hints' (selected), followed by an empty text field with browse and delete buttons.
 - Posterior File Name (optional):** An empty text field with browse and delete buttons.
 - ps2snp File (recommended):** A text field containing 'Axiom_PMRA.r1.ps2snp_map' with browse and delete buttons.

Using the Analysis Settings Fields

Follow the instructions below to create a new analysis configuration or edit a pre-populated field(s).

Sample QC Fields

1. Click the **Analysis File** drop-down button to select the appropriate XML file.
2. Click the **Prior Model File** Browse  button.
The Prior Model File window appears.
3. Navigate and select the appropriate file, then click **Open**.
Your newly assigned filename is displayed.
4. (Optional) Click the **SNP List File** Browse  button.
The SNP List File window appears.
5. Navigate and select the appropriate file, then click **Open**.
Your newly assigned filename is displayed.
6. (Optional) Click the **Gender File** Browse  button.
The Gender File window appears.
7. Navigate and select the appropriate file, then click **Open**.
Your assigned filename is displayed.
8. (Optional) Click the **Hints/Inbred File** Browse  button.
The Hints/Inbred File window appears.
9. Navigate and select the appropriate file, then click **Open**.
Your newly assigned path is displayed.
10. Click the either the **Inbred** or **Hints** radio button.

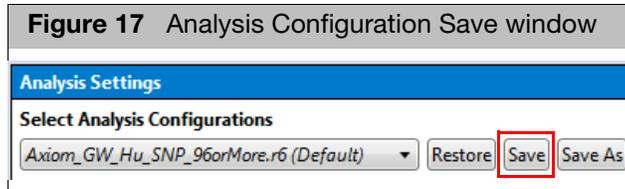
Genotyping Fields

1. Click the **Analysis File** drop-down button to select the appropriate XML file.
2. Click the **Prior Model File** Browse  button.
The Prior Model File window appears.
3. If multi-allelic probesets are available, the following fields appear:
 - Multi-allele Background Prior Model File
 - Multi-allele Pairwise Prior Model File
 - Multi-allele Prior Model File
 - If you want to change the currently displayed Model file, click the appropriate Multi-allele Browse  button to select a different Model file.
4. Navigate and select the appropriate file, then click **Open**.
Your newly assigned filename is displayed.
5. (Optional) Click the **SNP List File** Browse  button.
The SNP List File window appears.
6. Navigate and select the appropriate file, then click **Open**.
Your newly assigned filename is displayed.
7. (Optional) Click the **Gender File** Browse  button.
The Gender File window appears.

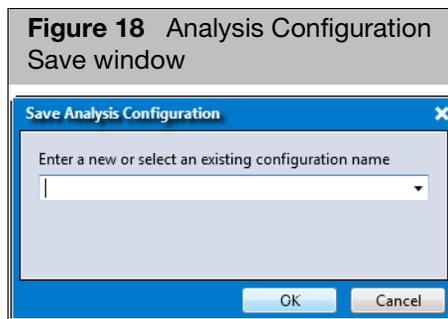
8. Navigate and select the appropriate file, then click **Open**.
Your assigned filename is displayed.
9. (Optional) Click the **Hints/Inbred File** Browse  button.
The Hints/Inbred File window appears.
10. Navigate and select the appropriate file, then click **Open**.
Your assigned filename is displayed.
11. Click either the **Inbred** or **Hints** radio button.
12. (Optional) Click the **Posterior File Name** Browse  button.
The Posterior File Name window appears.
13. Navigate to a location where you want to save your posterior file, enter a name, then click **Open**.
Your assigned filename is displayed.
14. Click the **ps2snp File** Browse  button.
The ps2snp File window appears.
15. Navigate to your ps2snp-file location, then click **Open**.
Your newly assigned filename is displayed.
16. (Optional) If the library package supports it, click the **Genotype Frequency** Browse button.
The Genotype Frequency window appears.
17. Navigate to a location for your genotype frequency file, enter a name, then click **Open**.
Your assigned filename is displayed.

Saving your Analysis Configuration

1. After editing your Analysis Configuration settings, click **Save** (top of Analysis Setting pane). (Figure 17)



If the Analysis Configuration that came with the library package already exists, then a **Save Analysis Configuration** window appears. (Figure 18)



2. Enter a different configuration name, then click **OK**.
Your new Analysis Configuration name is saved and available for use in the **Select Analysis Configuration** drop-down menu.

Modifying an Existing Analysis Configuration

1. Click the **Select Analysis Configuration** drop-down, then click to select the saved analysis configuration you want to modify.

Do one or more of the following to modify an existing analysis configuration:

- Click the applicable File field's **Browse** button to navigate to a different location, then click **Open** to reassign its path.
- If needed, click a File field's  button to delete a displayed path setting.
- Click the **Restore** button to return to the last saved values of the analysis configuration file.
- Click the **Save** button to overwrite your previously saved configuration
- Click the **Save As** button to save your modified configuration with a different name. [Recommended]

Setting Up Threshold Settings

The settings shown in the **Threshold Setting** pane (Figure 19) are based on the Mode (Workflow) you selected.

For Sample QC and SNP QC name definitions, see "Threshold Names" on page 149.

Figure 19 Automated QC Mode Threshold Settings pane example

Threshold Settings

Select Threshold Configurations
Human (Default) Restore Save Save As

Sample QC

Name	Settings
DQC	≥ 0.82
QC call_rate	≥ 97
Percent of passing sampl...	≥ 95
Average call rate for pass...	≥ 98.5

SNP QC

Name	Settings
species-type	Human
cr-cutoff	≥ 95
fld-cutoff	≥ 3.6
het-so-cutoff	≥ -0.1
het-so-otv-cutoff	≥ -0.3
hom-ro-1-cutoff	≥ 0.6
hom-ro-2-cutoff	≥ 0.3
hom-ro-3-cutoff	≥ -0.9
hom-ro	true
hom-het	true
num-minor-allele-cutoff	≥ 2
priority-order	Change List Order PolyHighResolution, NoMi...
recommended	Checklist PolyHighResolution, NoMinorHom,...

Customizing Thresholds

Click the **Select Threshold Configuration** drop-down (Figure 20) to select an appropriate Default Threshold for your starting point. Note: The comparison signs/operators are preset and cannot be changed.

Figure 20 Select Threshold Configuration

Select Threshold Configurations

- Diploid (Default)
- Diploid (Default)
- Human (Default)
- Polyploid (Default)

Sample QC

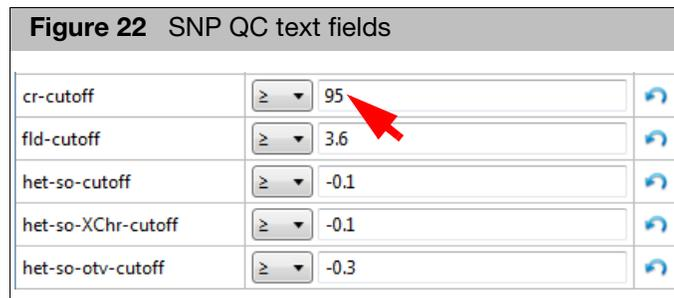
All the Sample QC Threshold Settings are populated with default values.

1. Click inside each text field to enter a different value, as shown in [Figure 21](#).
Click the text field's  button to return its value back to its last saved value within the threshold configuration file.



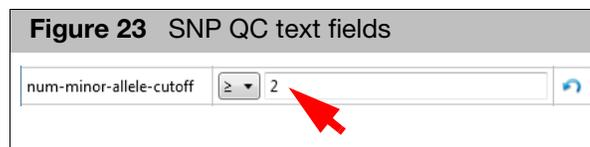
SNP QC

1. Click the **species-type** drop-down menu to select a different species type.
2. Click inside each text field to enter a different value, as shown in [Figure 22](#). Note: General Rule: The het-so-otv-cutoff should be less or equal to het-so-cutoff and het-so-XChr-cutoff.



Note: To return a value back to its last saved value, click the text field's  button.

3. Use the **hom-ro** and **hom-het** drop-down menus to change their True or False values.
4. Click inside the **num-minor-allele-cutoff** text field to enter a different value, as shown in [Figure 23](#).



5. The priority-order option enables you to change the order of categories when determining which probesets are selected as the best probeset for a SNP. To change the priority-order of your SNP QC Metric, click [Change List Order](#).

The following window appears: (Figure 24)

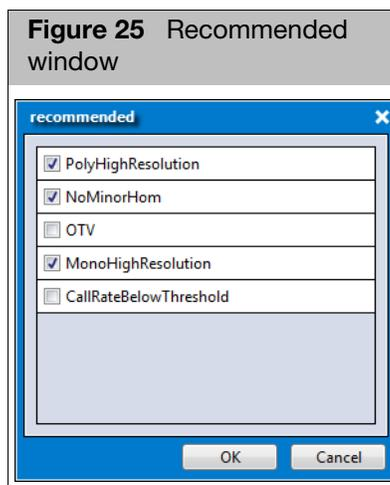


- Click and hold onto the selection you want to move, then drag and drop it into its new position. After you get the order of priority you want, click **OK**.

Note: To return the list back to its default priority, click the priority-order field's  button.

- Use the **recommended** checklist to choose the PS_Classification conversion types for your analysis. To change the recommended options, click **Checklist**.

The recommended window appears. (Figure 25)



- Click to check/uncheck the available recommended options, then click **OK**.

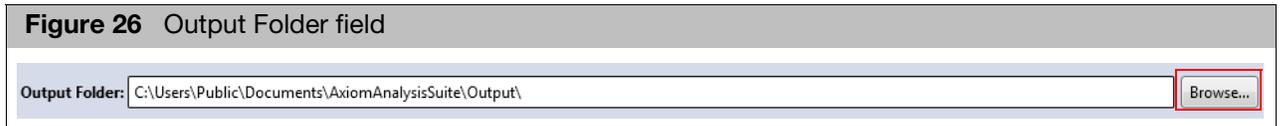
Note: If all recommended options are unchecked, the software uses the following default values:

- Human:** PolyHighResolution, NoMinorHom, MonoHighResolution
- Diploid:** PolyHighResolution, NoMinorHom, MonoHighResolution
- Polyploid:** PolyHighResolution

Assigning an Output Folder Path

Assigning a New Output Folder Path

1. Click the **Output Folder** path's **Browse** button. (Figure 26)



An Explorer window appears.

2. Navigate to the recommended path
C:\Users\Public\Documents\AxiomAnalysisSuite\Output, then click **Select Folder**.

Your selected output folder path is now displayed.

Adding Sub-Folders

Note: To better organize your output results, you can add sub-folders to your newly assigned output result path's folder.

To add sub-folders to your newly assigned result path's folder:

1. Click the Output Folder's **Browse** button to return to your assigned output path and/or folder.
2. In the Explorer window, click **New Folder**.
3. Enter a sub-folder name.
4. Click **Select Folder**.

The newly created sub-folder now appears in the output result information window.

5. Repeat the above steps 1-4 to add more sub-folders, then click **Select Folder**.

Assigning a Batch Name

The batch file is produced while your analysis is running and includes all the necessary files needed to view your analysis in the Axiom Analysis Suite Viewer.

1. Enter a name in the **Batch Name** field. (Figure 27)

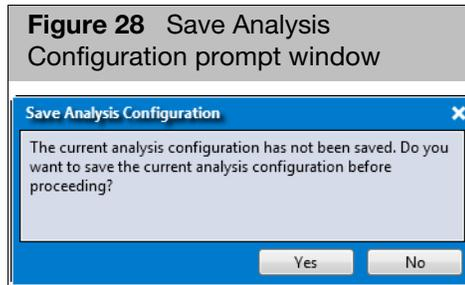
✗ IMPORTANT! Each Batch Name you create must be unique for the set of batches listed in the Dashboard window tab and unique within the same destination folder.



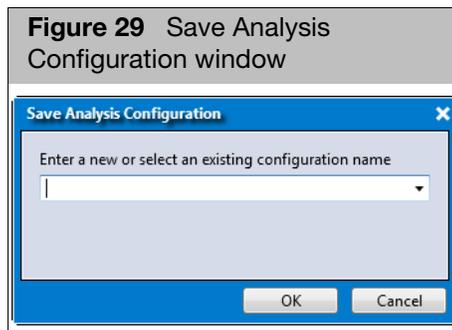
Note: A folder (with the same name as your entered batch name) is auto-generated during the analysis process. This folder includes all the necessary files needed to view your analysis results in the Viewer.

Running your Analysis

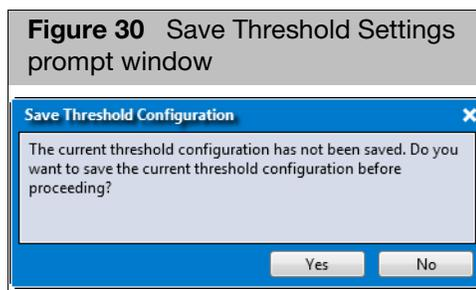
1. Click **Run Analysis**.
 - If you have not saved any changes to your configured Analysis Settings, a Save Analysis Configuration window appears. (Figure 28) Click **Yes**.



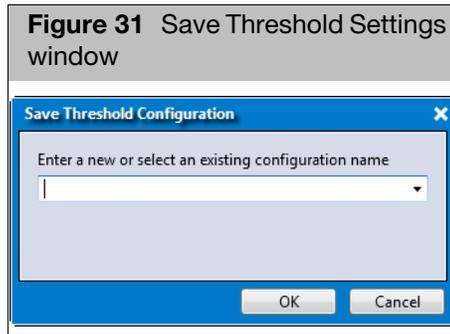
- Enter a new analysis name or use the drop-down to select a previously saved name, then click **OK**. (Figure 29)



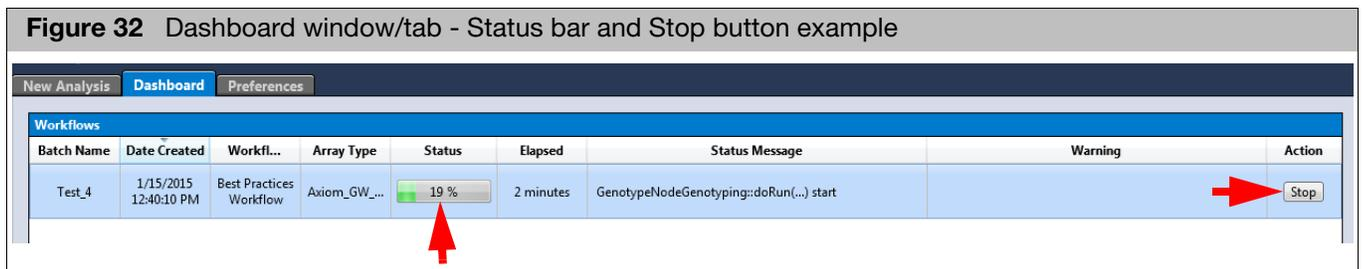
- If you have not saved any changes to your configured Threshold Settings, a Save Threshold Configuration window appears. (Figure 30) Click **Yes**.



- Enter a new threshold name or use the drop-down to select a previously saved name, then click **OK**.

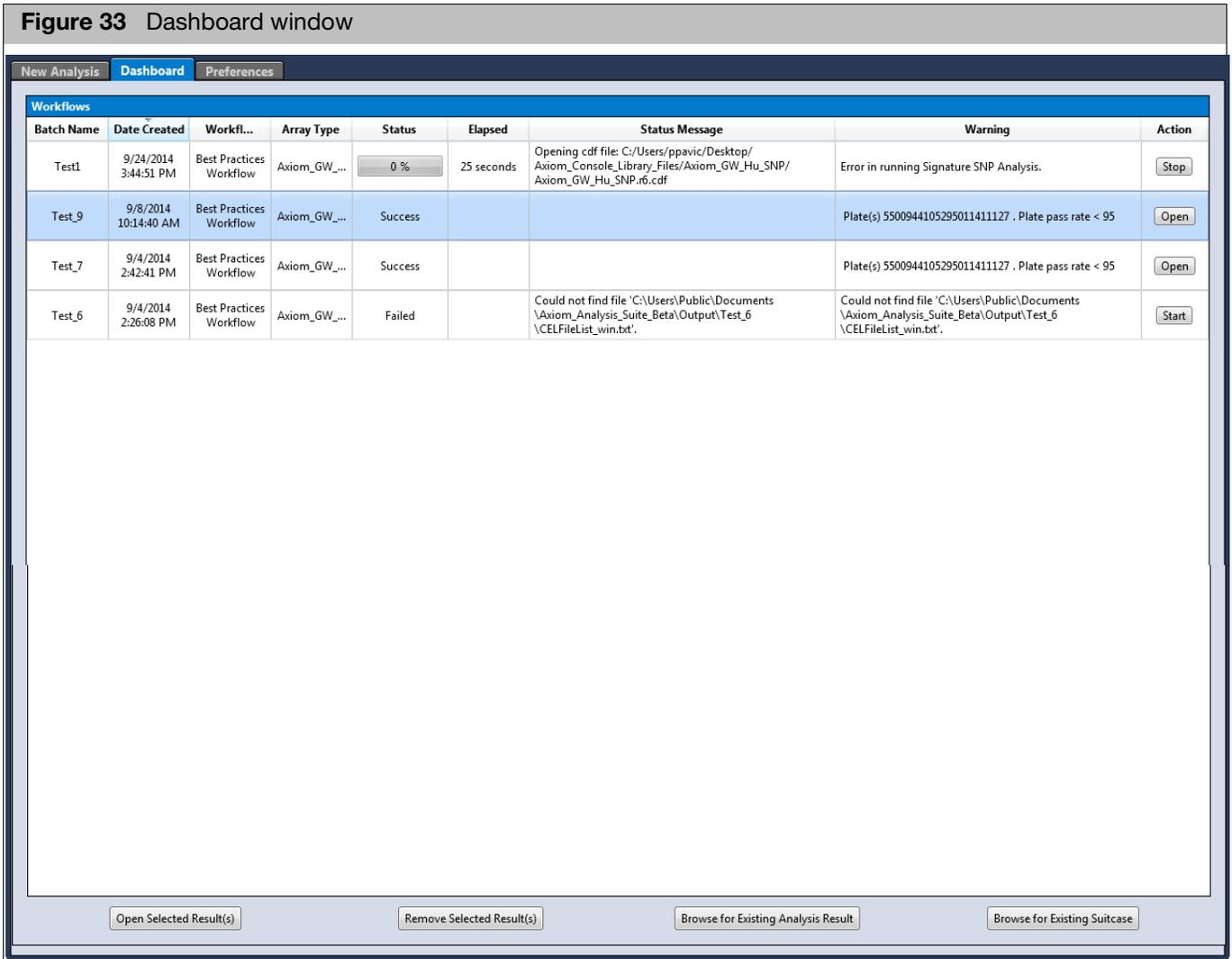


The Dashboard window/tab appears and shows the status of your running analysis. (Figure 32) Click **Stop** to cancel an analysis in progress.



Using the Dashboard Window Tab

The Dashboard tab window displays existing results. (Figure 33)



Open Selected Result(s)

Do one of the following to open a selected result:

- Click on the Action column's **Open** button to open a study.

Note: The **Open** button is relabeled **Start** if the Dashboard's Status column displays Failed. A workflow can fail if there is a processing error (for example, needed files are unavailable) or if no samples pass QC. If the Warning column indicates that no samples passed QC, yet you still want to open the study to review its QC results, use one of the alternate methods below to open a selected study.

- Single-click on a study, then click **Open Selected Result(s)**.
- Double-click on a study to open it.
- Right-click on a study, then click **Open**. Note: Click on any of the Workflow header columns to sort your listed results.

After a few moments, the Viewer opens and displays your study (as you last left it).

Remove Selected Result(s)

Do one of the following to remove a selected result:

- Single-click to highlight the analysis you want to remove, then click **Remove Selected Result(s)**.
- Right-click on the highlighted analysis, then click **Remove from List**.

Viewing the Results Folder in Windows Explorer

1. Right-click on the highlighted analysis, then click **Open in Windows Explorer**. Your Analysis Results folder now appears in the Explorer window.

Browsing For Existing Analysis Results

The Axiom Analysis Suite Dashboard displays/stores your results for each user profile. Note: Navigate to the Workflows sub-folder to locate other User Profile folders (that may contain more recent results).

Do the following if a study is not listed on the Dashboard:

1. Click **Browse for Existing Analysis Result**.
A Select Analysis Result File window appears.
2. Click on a recent analysis, then click **Select Folder**.
After a few moments, your analysis opens as you last left it.
3. After your analysis has successfully completed. click **Open** (Figure 34).
Alternatively, click to highlight the completed analysis, then click **Open Selected Result(s)**.



The Axiom Analysis Suite Viewer appears.

For instructions on how to use the Viewer, continue to [Chapter 3](#).

Browsing for Existing Suitcases

In AxAS v1.0, a suitcase file was auto-generated during an analysis and included all the necessary files needed to view your analysis in the Viewer.

For AxAs v1.1 and newer, the suitcase has been replaced with a higher performing batch name folder. Your suitcase file must be converted to this new folder format before it can be opened in the newly installed Viewer.

Do the following to convert your suitcase file to a batch name folder:

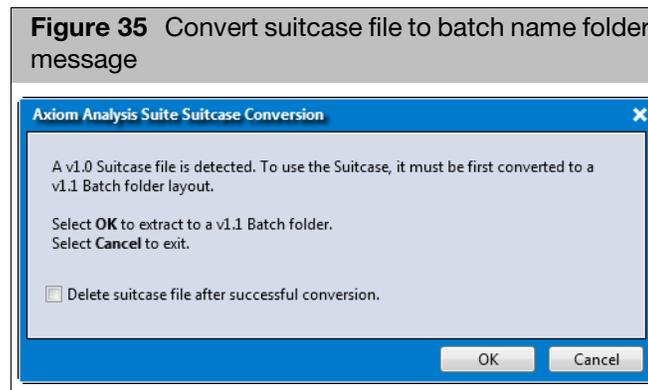
1. Click **Browse for Existing Suitcase**.

A Select Analysis Result File window appears.

2. Click to highlight a suitcase file, then click **Open**.

An **Axiom Analysis Suite Suitcase Conversion** message window appears.

(Figure 35)



3. If you want to retain your v1.0 suitcase file for archiving purposes, leave the **Delete suitcase file after successful conversion** check box unchecked. Click on this check box if you want your suitcase file to be auto-deleted after it is converted.

4. Click **OK**.

Allow a few moments for your suitcase file to convert to the v1.1 batch name folder format.

The Axiom Analysis Suite Viewer appears.

For instructions on how to use the Viewer, continue to [Chapter 3](#).

Opening an Analysis from Windows Explorer

You can open an existing analysis directly from Windows Explorer, instead of through the application's Dashboard window tab. To do this:

1. Right-click on an analysis batch folder and select **Open in Axiom Analysis Suite Viewer**.

A Select Profile window appears.

2. At the Select Profile window, enter a new or select an existing profile name.

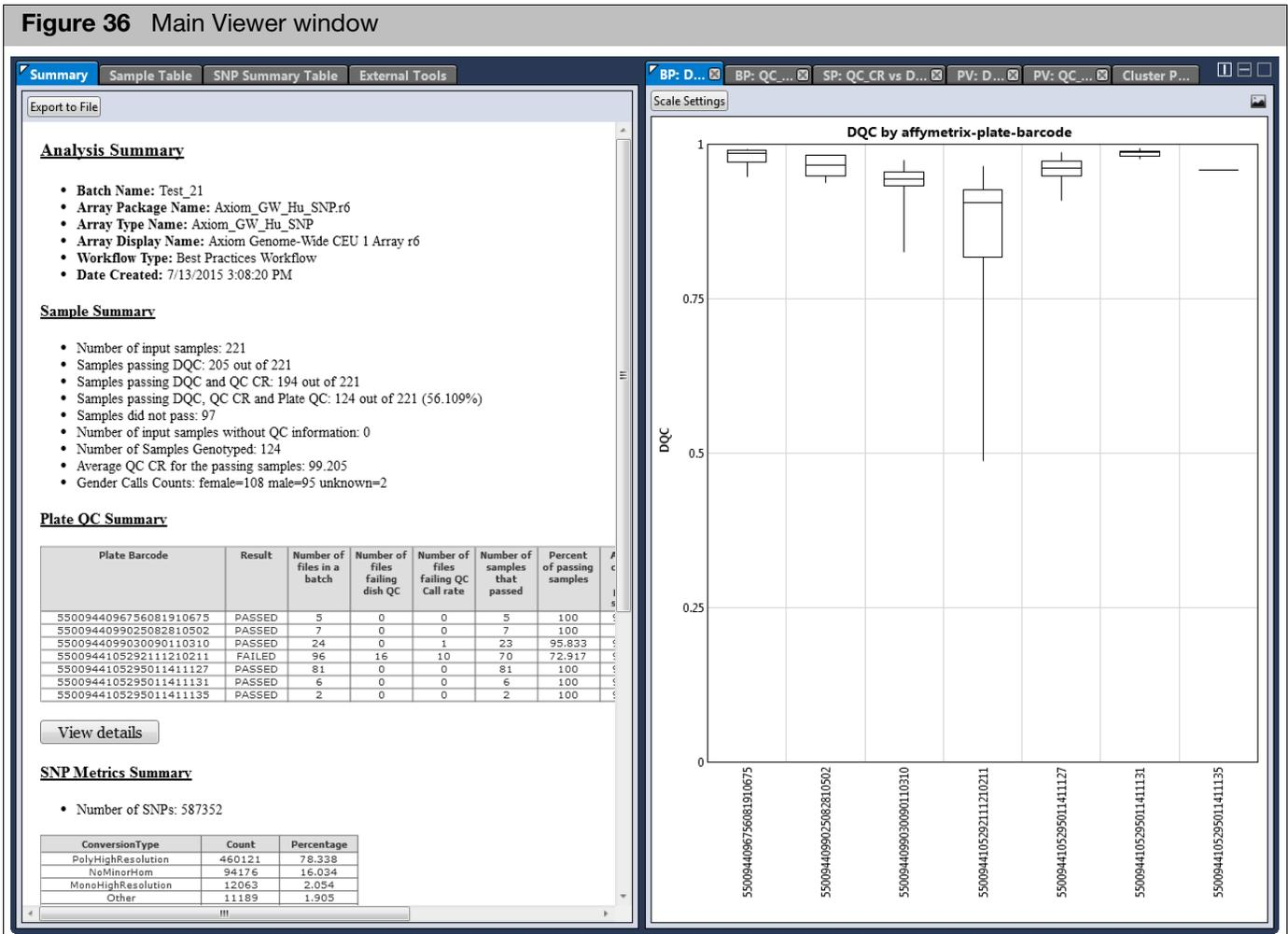
If your selected folder is a valid analysis folder, the application's Viewer opens as it normally would.

3

The Viewer: Summary Window and Sample Table

After setting up and successfully running an analysis, as described in [Chapter 2](#), the Axiom Analysis Suite Viewer opens. (Figure 36)

Figure 36 Main Viewer window



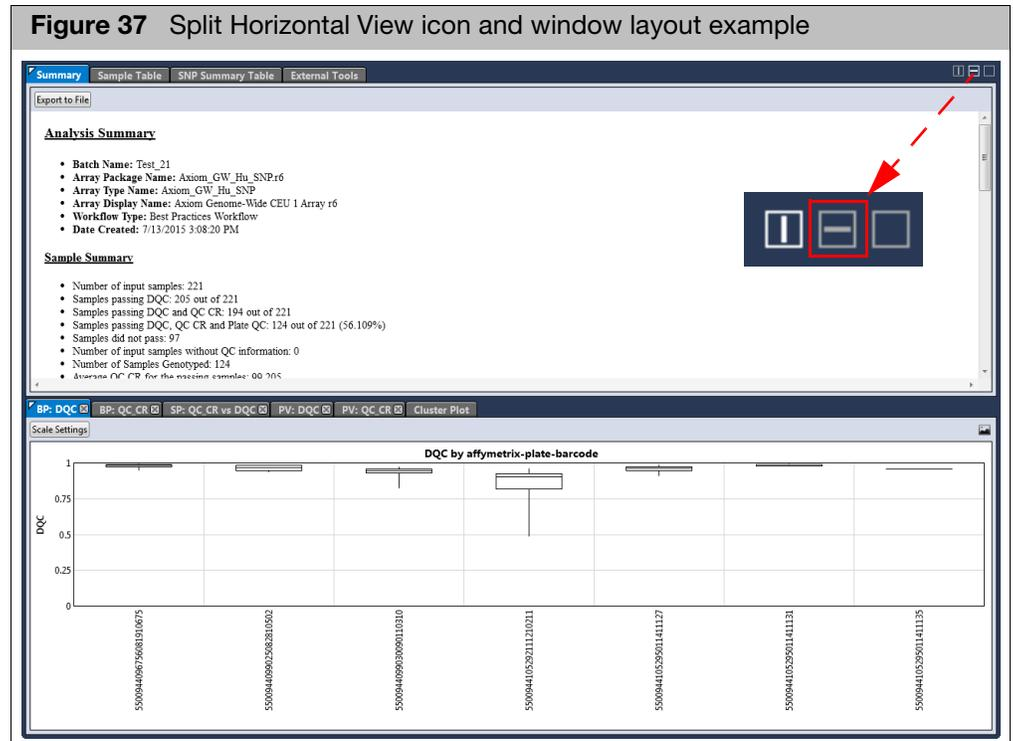
Viewing Options

As shown in [Figure 36](#) on page 38, the Viewer (by default) displays a **side-by-side** split-screen configuration.

Split-Screen Options

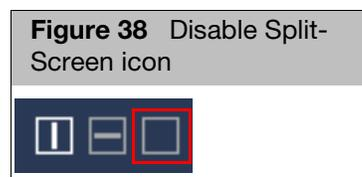
To change side by side split-screen to a top and bottom configuration:

1. Click the **Horizontal Split** icon. ([Figure 37](#))

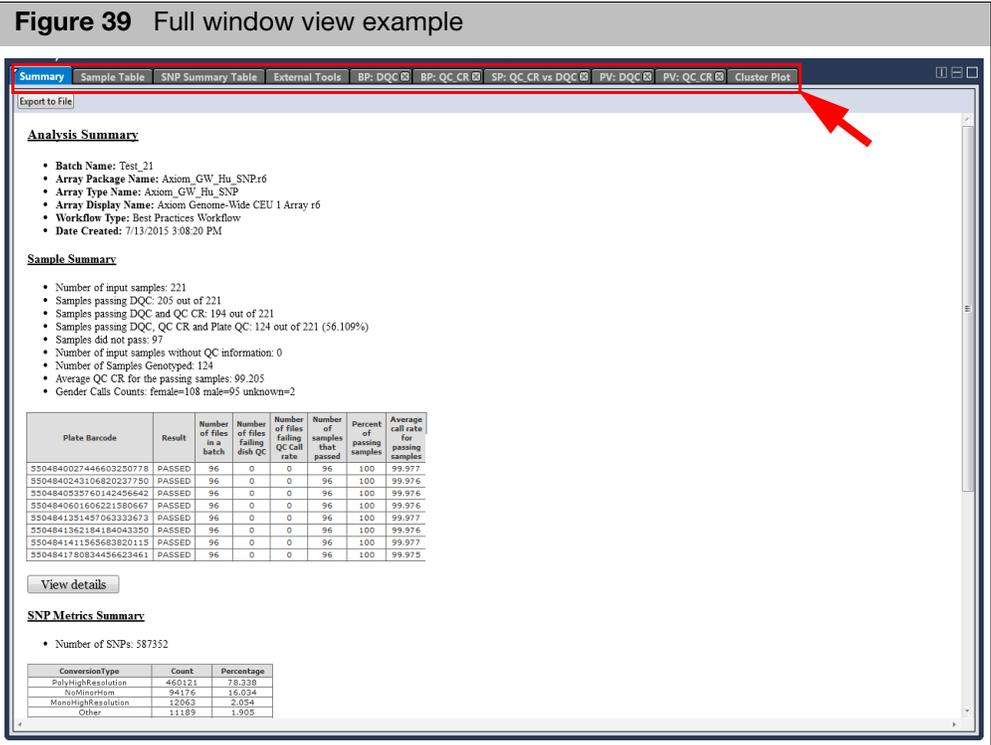


To disable the split-screen:

1. Click the **Disable Split-Screen** icon. ([Figure 38](#))



The split-screen becomes 1 window. ([Figure 39](#))



2. Click on any window tab (Figure 39) to view it in full window mode.

To return to the default side by side split-screen configuration:

1. Click the Vertical Split icon. (Figure 40)

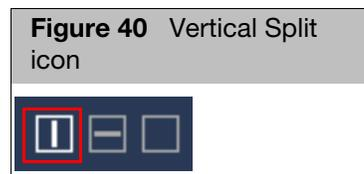
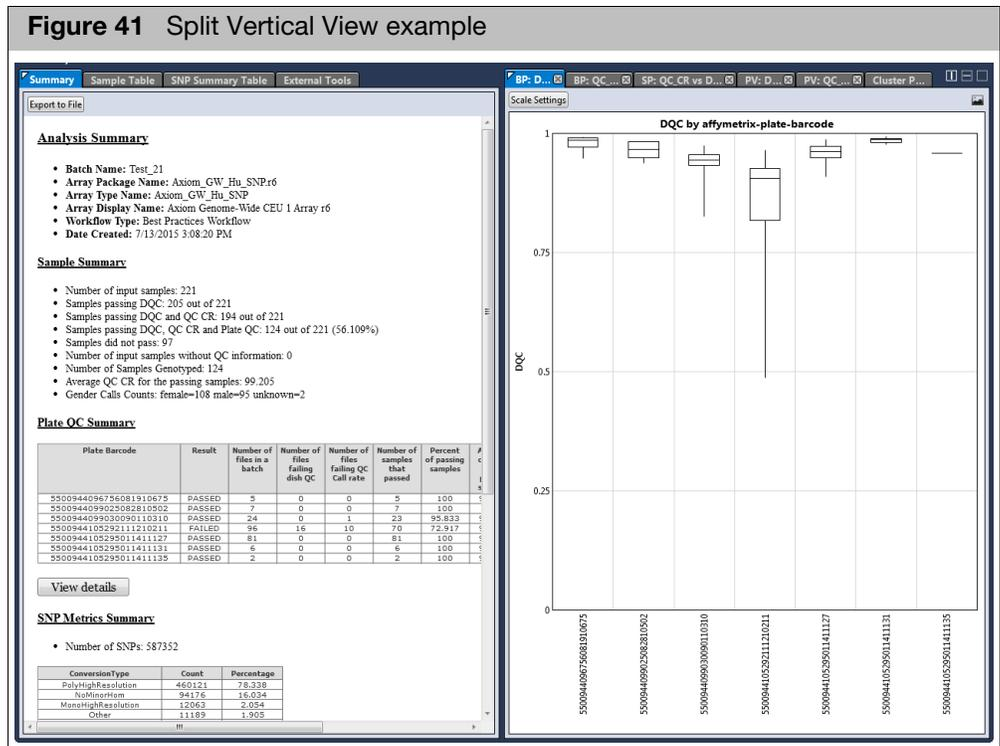


Figure 41 Split Vertical View example



Changing a Tab Window to a Full Screen Windows

To toggle a tab window to full screen:

1. Locate the tab you want to make full screen.
2. Click on a tab's white triangle graphic. (Figure 42)



The window tab is now a window.

3. Double-click anywhere along the top of the window to change it to full screen.

To toggle a full screen window to its default tab window:

1. Double-click anywhere along the top of the window.
2. Click the X button (top right) to close the window.

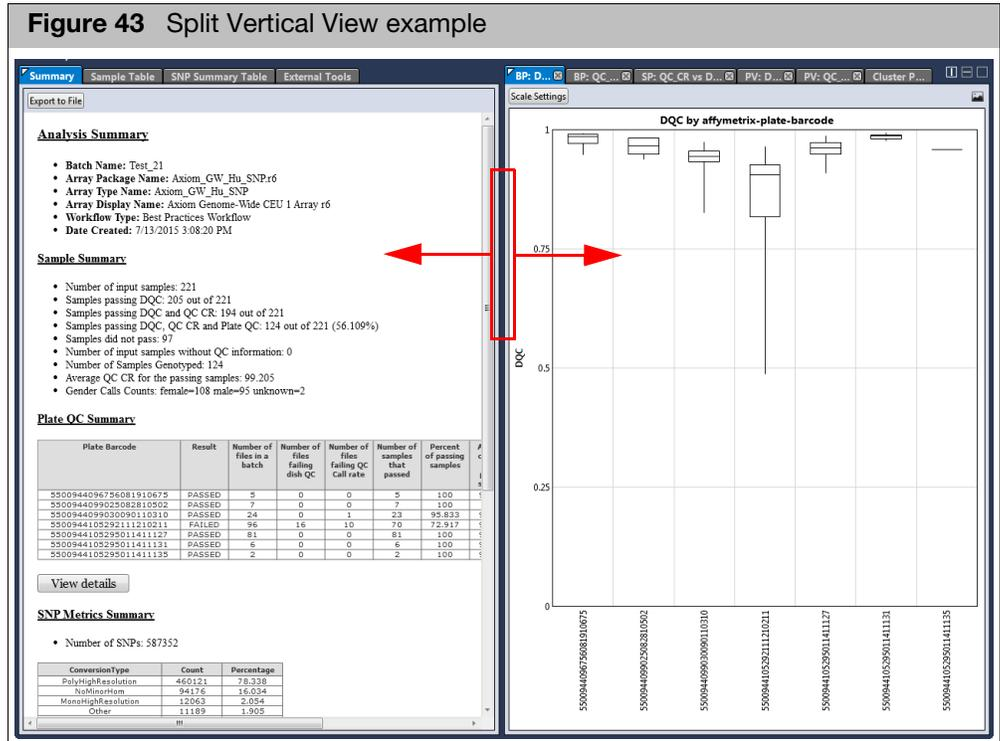
The window returns to its default tab window and position.

Adjusting the Window Size

To change the size of a window pane:

1. Click, hold, then drag the edge of the window pane (Figure 43) to resize it.

Figure 43 Split Vertical View example



Summary Window/Tab

The Summary window/tab (Figure 44) displays a summary snapshot of your analysis, including detailed threshold values, and tables based on your analysis.

Data Analysis Summary

Note: Each workflow type reports different information within the Analysis Summary window. Figure 44 is an example of a Best Practices workflow.

Figure 44 Summary window tab

Export to File

Analysis Summary

- **Batch Name:** Test_21
- **Array Package Name:** Axiom_GW_Hu_SNP:r6
- **Array Type Name:** Axiom_GW_Hu_SNP
- **Array Display Name:** Axiom Genome-Wide CEU 1 Array r6
- **Workflow Type:** Best Practices Workflow
- **Date Created:** 7/13/2015 3:08:20 PM

Sample Summary

- Number of input samples: 221
- Samples passing DQC: 205 out of 221
- Samples passing DQC and QC CR: 194 out of 221
- Samples passing DQC, QC CR and Plate QC: 124 out of 221 (56.109%)
- Samples did not pass: 97
- Number of input samples without QC information: 0
- Number of Samples Genotyped: 124
- Average QC CR for the passing samples: 99.205
- Gender Calls Counts: female=108 male=95 unknown=2

Plate QC Summary

Plate Barcode	Result	Number of files in a batch	Number of files failing dish QC	Number of files failing QC Call rate	Number of samples that passed	Percent of passing samples
5500944096756081910675	PASSED	5	0	0	5	100
5500944099025082810502	PASSED	7	0	0	7	100
5500944099030090110310	PASSED	24	0	1	23	95.833
5500944105292111210211	FAILED	96	16	10	70	72.917
5500944105295011411127	PASSED	81	0	0	81	100
5500944105295011411131	PASSED	6	0	0	6	100
5500944105295011411135	PASSED	2	0	0	2	100

View details

SNP Metrics Summary

- Number of SNPs: 587352

ConversionType	Count	Percentage
PolyHighResolution	460121	78.338
NoMinorHom	94176	16.034
MonoHighResolution	12063	2.054
Other	11189	1.905
CallRateBelowThreshold	7568	1.288
Hemizygous	1971	0.336
OTV	264	0.045

Sample QC Thresholds

- DQC: ≥ 0.81
- QC call_rate: ≥ 97
- Percent of passing samples: ≥ 88
- Average call rate for passing samples: ≥ 98.5

SNP QC Thresholds

- species-type: Diploid
- cr-cutoff: ≥ 94
- fld-cutoff: ≥ 3.6
- het-so-cutoff: ≥ -0.1

Export to File: Click this button to export the Summary report as a PDF file.

Analysis Summary: Contains information about the array type, the workflow run and the date processed

Sample Summary: Breaks down the sample QC for your analysis run and displays the number that pass each of your QC Thresholds. In addition, it provides the average QC Call Rate (CR) and breakdown of the genders found within your batch of samples.

Plate QC Summary: Contains sample QC information for each plate including the number samples failing DQC, QC Call Rate, the Percent of passing samples, and the average Call Rate for your passing samples.

SNP Metrics Summary: This section contains a summary of the categorization of the SNPs in the analysis by PS_Classification. For more information on these categories see *“Regenerate SNP Metrics” on page 82.*

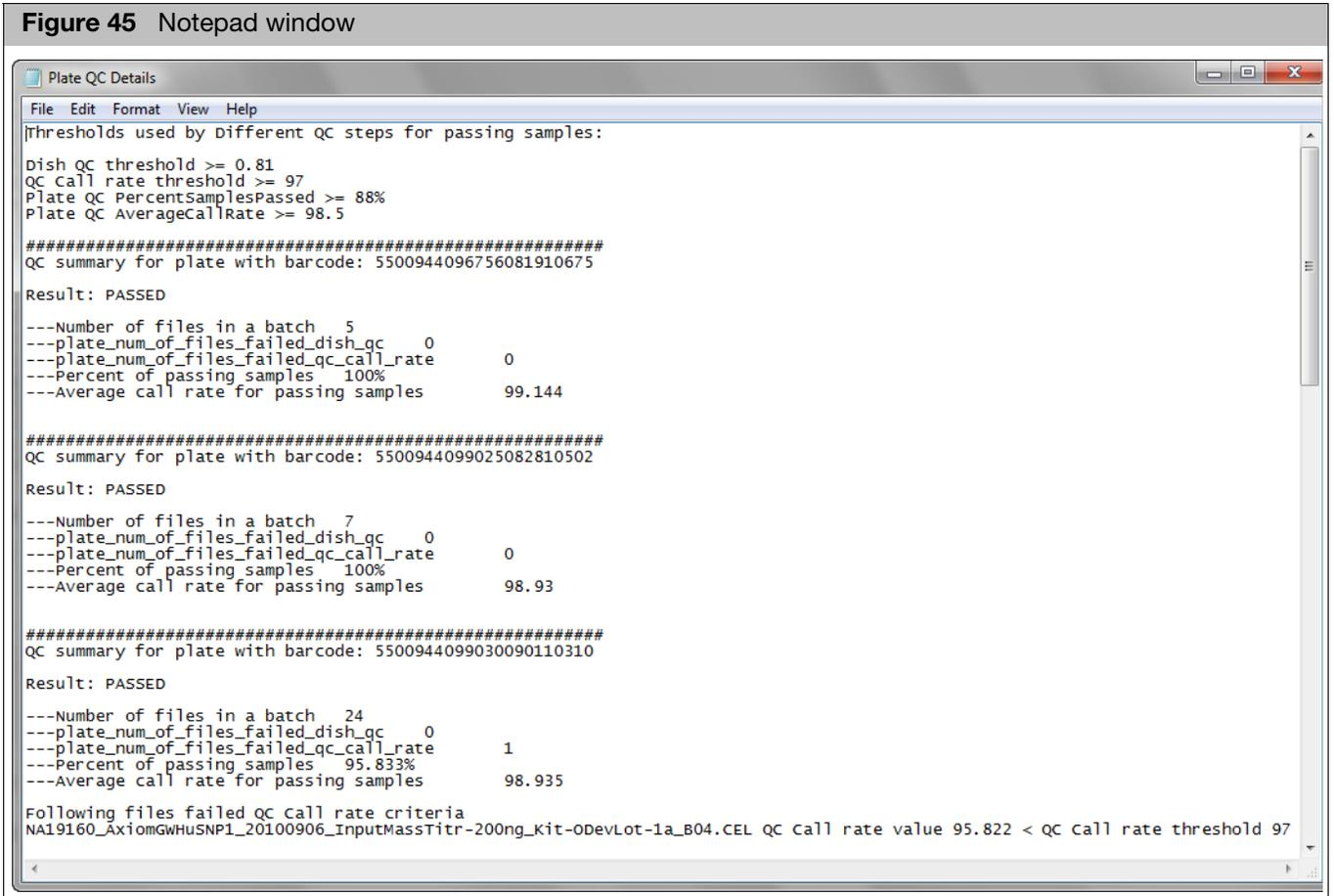
Sample QC Thresholds: Displays the Sample QC Thresholds used for your analysis run and their associated SNP QC Metrics.

SNP QC Thresholds: Displays the Thresholds used for your analysis run and their associated SNP QC Metrics.

Viewing the Plate Barcode Table Details

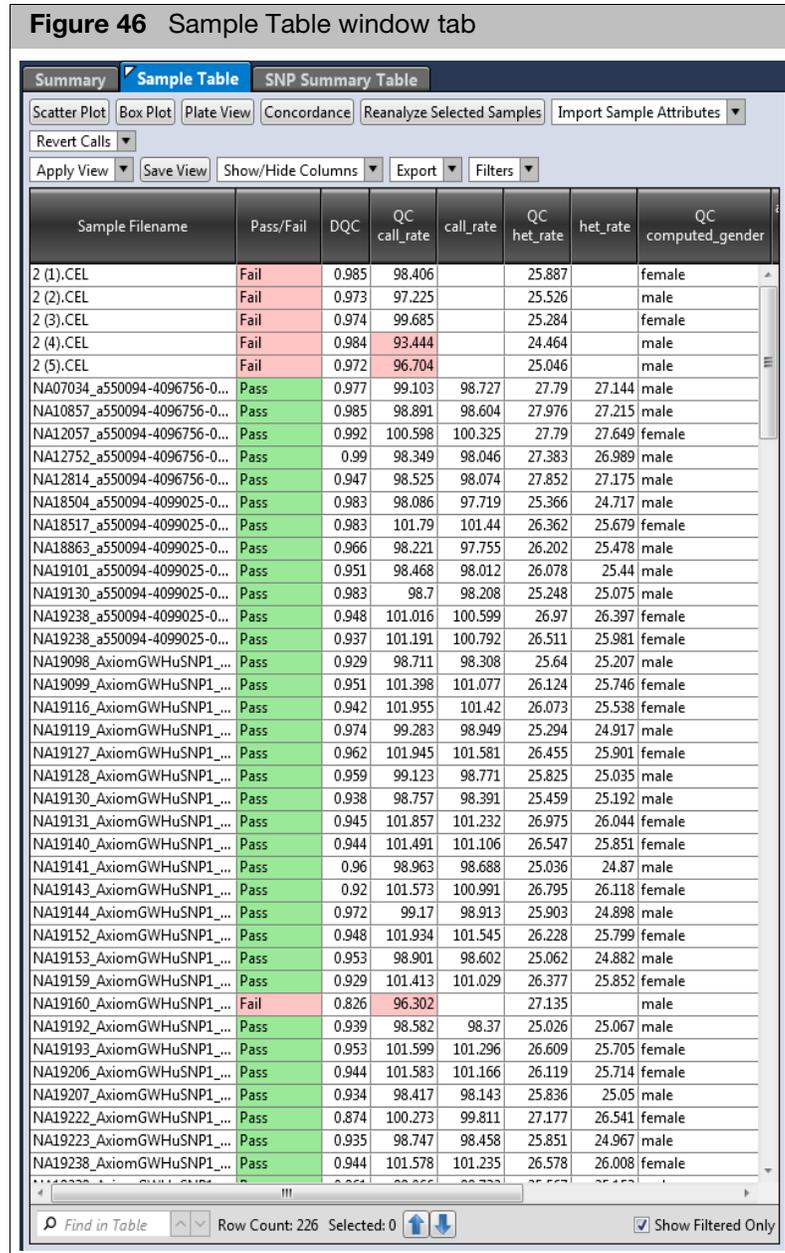
1. In the Summary window tab, click **View Details**. (Figure 44)
A window opens and displays a text file version of your Sample QC information (by plate). (Figure 45)

Figure 45 Notepad window



Sample Table

Note: Depending on the Threshold values you set (prior to running your analysis), color-coded Pass or Fail cells may appear in the table, as shown in Figure 46.



Importing Sample Attributes

To import sample attributes into your Sample Table:

1. Click the **Import Sample Attributes** drop-down.
2. Click to select either **Import from ARR Files** or **Import from CSV/Tab-Delimited Text File**.

An Explorer window appears.

✗ IMPORTANT! Your text-based CEL file must start with the header *Sample Filename* and include the full CEL file name, as shown in [Figure 47](#).

Figure 47 Tab-delimited text CEL file example shown in Excel

	A	B	C
1	Sample Filename	String Name	Number Name
2	NA18961_AxiomGWHuSNP1_20101007_TrainingMix_Rtstab-DL1a-Omo_E09.CEL	A1	1
3	NA12760_AxiomGWHuSNP1_20101007_TrainingMix_Rtstab-DL1a-Omo_B06.CEL	A2	1
4	NA19160_AxiomGWHuSNP1_20100906_InputMassTitr-200ng_Kit-ODevLot-1a_B04.CEL	A3	1
5	NA18960_AxiomGWHuSNP1_20101007_TrainingMix_Rtstab-DL1a-Omo_E08.CEL	A4	1
6	NA12814_AxiomGWHuSNP1_20101007_TrainingMix_Rtstab-DL1a-Omo_B09.CEL	A5	1
7	NA19210_AxiomGWHuSNP1_20101007_TrainingMix_Rtstab-DL1a-Omo_F12.CEL	A6	1

3. Navigate to the applicable file location, then click **Open**.

Column Headers

The default Sample Table column view is as shown. ([Figure 48](#))

Figure 48 Default Sample Table Columns

Sample Filename	Pass/Fail	DQC	QC call_rate	call_rate	QC het_rate	het_rate	QC computed_gender	affymetrix-plate-barcode	affymetrix-plate-peg-wellposition	Average call rate for passi...	Percent of passing...
-----------------	-----------	-----	--------------	-----------	-------------	----------	--------------------	--------------------------	-----------------------------------	--------------------------------	-----------------------

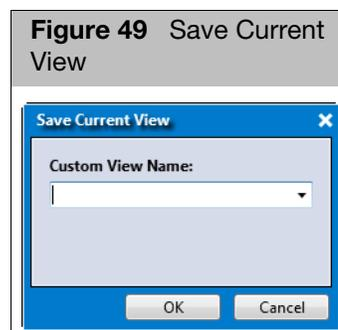
To show or hide table columns:

1. Click the **Show/Hide Columns** drop-down menu.
2. Click each available column name's check box to show it or remove it from the table. See "[Annotations and Columns](#)" on page 159 for their definitions.
3. Click outside the **Show/Hide Columns** drop-down menu to close it.

To save your customized Sample Table column view:

1. Click **Save View**.

The Save Current View window appears. ([Figure 49](#))



2. Enter a name for your custom table view, then click **OK**.

Your newly saved name is now added to the **Apply View** drop-down menu.

To show ALL available columns within the Sample Table:

1. Click the **Apply View** drop-down menu, then select **All Columns View**.

Rearranging Columns

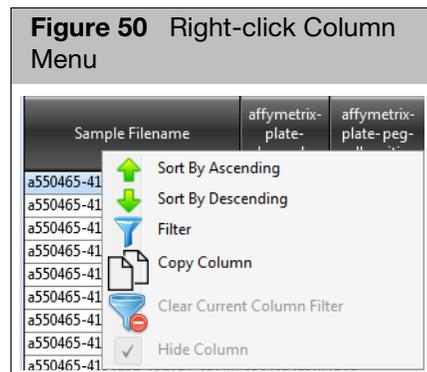
1. Click on a column you want to move.
2. Drag it (left or right) to its new location.
3. Release the mouse button.

The column is now in its new position.

Sorting Columns

1. Select a column, then right-click on it.

A right-click menu appears. (Figure 50)



2. Click to select either **Sort By Ascending** (A-Z) or **Sort By Descending** (Z-A).

Single-Click Sorting Method

1. Single-click on a column header to sort its data in an ascending order. Single-click on the same column header to sort its data in a descending order

Hiding the Column

1. Select the column you want to hide from the table, then right-click on it.

A right-click menu appears. (Figure 50)

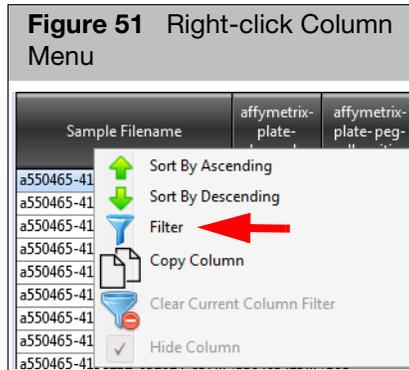
2. Click the **Hide Column** check box to remove it from the table.

Filtering Column Data

Note: All Sample Table columns are filterable.

Adding Filters (Method 1)

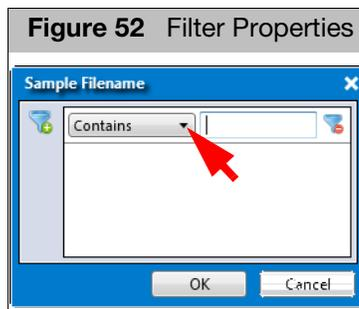
1. Select a column, then right-click on it.
The following window appears: (Figure 51)



2. Click **Filter**.

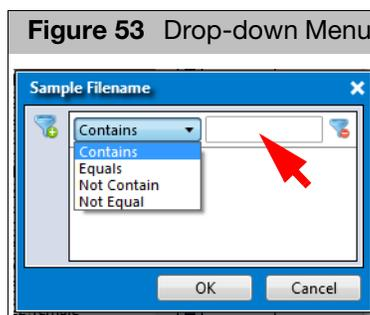
Text-based Columns

If the column you want to filter contains text-based data, the Contains drop-down menu appears. (Figure 52)

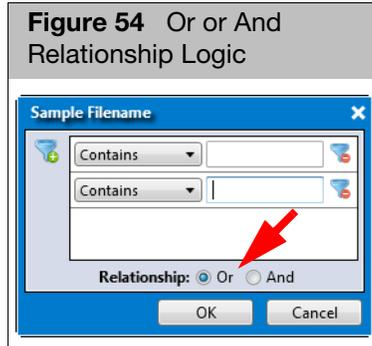


To apply a filter to a text-based column:

1. Click the **Contains** drop-down menu to select a filtering property. (Figure 53)



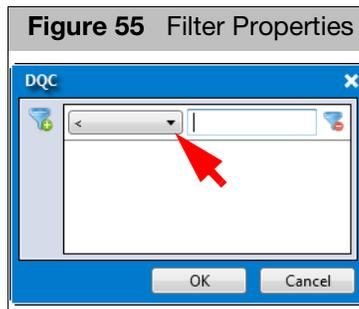
2. Click inside the text entry box to enter a value. (Figure 53)
3. OPTIONAL: Click  to add additional filters.



4. Click the **Or** or **And** radio button to choose **Or** or **AND** relationship logic. (Figure 54)
5. Repeat steps 1-4 as needed.
6. To remove a filter(s), click .

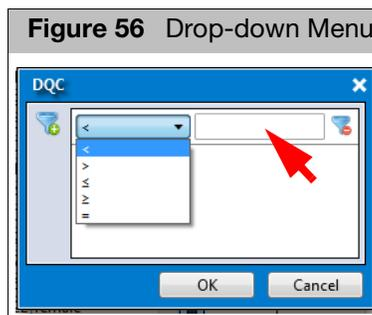
Numeric Data Columns

If the column you want to filter contains numeric data, a symbol drop-down menu appears. (Figure 55)



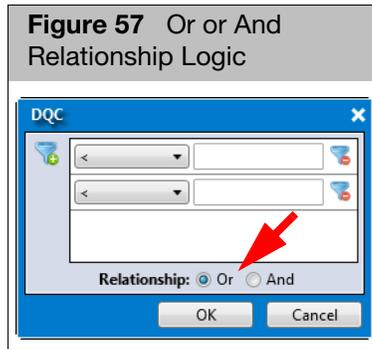
To apply a filter to a value-based column:

1. Click the **Symbol Value** drop-down menu to select the filtering symbol you want. (Figure 56)



2. Click inside the text entry box to enter the value(s). (Figure 56)

3. OPTIONAL: Click  to add filter(s).



4. Click the **Or** or **And** radio button to choose **Or** or **AND** relationship logic. (Figure 57)
5. If needed, repeat steps 1-4.
6. Click **OK**.
To remove a filter(s), click .

Showing Filtered Data Only

- Click the **Show Filtered Only** check box Show Filtered Only to show only the data that passes the filters.

Uncheck this box to show all data, including data that did not pass your filter criteria setting(s). In this mode, data that passes the filter appears in light gray, as shown in [Figure 58 on page 51](#).

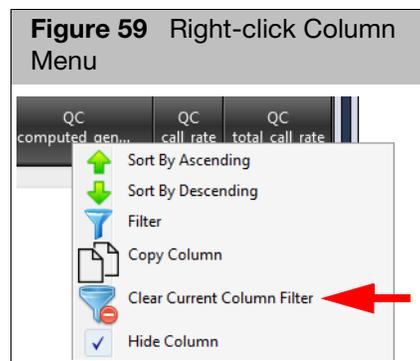
Figure 58 Sample Table window tab - Show Filter Only unchecked example

Sample Filename	Pass/Fail	DQC	QC call_rate	call_r...	QC het_rate	het_r...	QC computed_gen...
NAI1840_AxiomGWHuSNP1_...	Pass	0.871	100.371	99.964	29.57	29.226	female
NAI2875_AxiomGWHuSNP1_...	Pass	0.833	99.479	98.943	30.292	29.594	female
NAI8577_AxiomGWHuSNP1_...	Pass	0.847	99.381	99.02	26.831	26.318	female
NAI8593_AxiomGWHuSNP1_...	Pass	0.902	99.665	99.318	26.403	25.716	female
NAI8942_AxiomGWHuSNP1_...	Pass	0.869	100.289	99.987	26.599	25.908	female
NAI8949_AxiomGWHuSNP1_...	Pass	0.925	99.175	98.832	25.913	24.988	female
NAI9099_AxiomGWHuSNP1_...	Pass	0.851	100.397	99.973	27.079	26.676	female
NAI9222_AxiomGWHuSNP1_...	Pass	0.874	100.273	99.811	27.177	26.541	female
2877_AxiomGWHuSNP1_2010...	Pass	0.982	102.424	102.115	26.847	25.092	female
2878_AxiomGWHuSNP1_2010...	Pass	0.972	102.357	101.935	25.49	24.78	female
2881_AxiomGWHuSNP1_2010...	Pass	0.951	101.96	101.588	28.585	27.786	female
2884_AxiomGWHuSNP1_2010...	Pass	0.955	102.089	101.636	26.083	25.184	female
2885_AxiomGWHuSNP1_2010...	Pass	0.969	102.213	101.81	27.749	27.9	female
2888_AxiomGWHuSNP1_2010...	Pass	0.945	101.676	101.279	28.471	28.032	female
2889_AxiomGWHuSNP1_2010...	Pass	0.986	102.563	102.292	28.678	27.89	female
2890_AxiomGWHuSNP1_2010...	Pass	0.97	102.352	101.956	26.393	25.224	female
2893_AxiomGWHuSNP1_2010...	Pass	0.965	102.357	101.93	28.781	28.135	female
2894_AxiomGWHuSNP1_2010...	Pass	0.967	102.233	101.855	25.897	25.244	female
2895_AxiomGWHuSNP1_2010...	Pass	0.978	102.527	102.255	25.676	24.955	female
2897_AxiomGWHuSNP1_2010...	Pass	0.965	102.553	102.1	26.248	25.142	female
2899_AxiomGWHuSNP1_2010...	Pass	0.969	102.373	101.972	28.796	28.016	female
2900_AxiomGWHuSNP1_2010...	Pass	0.953	102.094	101.667	28.177	28.08	female
2901_AxiomGWHuSNP1_2010...	Pass	0.979	102.357	101.977	28.409	27.583	female
2902_AxiomGWHuSNP1_2010...	Pass	0.979	102.486	102.163	26.955	25.39	female
2903_AxiomGWHuSNP1_2010...	Pass	0.975	102.507	102.082	28.961	28.181	female
2905_AxiomGWHuSNP1_2010...	Pass	0.965	102.285	101.844	28.09	27.9	female
2907_AxiomGWHuSNP1_2010...	Pass	0.95	102.104	101.603	26.119	25.442	female
2908_AxiomGWHuSNP1_2010...	Pass	0.951	102.146	101.701	28.93	27.792	female
2909_AxiomGWHuSNP1_2010...	Pass	0.963	102.383	102.008	25.418	24.772	female
2915_AxiomGWHuSNP1_2010...	Pass	0.961	102.151	101.827	28.817	28.035	female
2916_AxiomGWHuSNP1_2010...	Pass	0.985	102.481	102.216	26.063	24.87	female
2917_AxiomGWHuSNP1_2010...	Pass	0.956	102.12	101.717	28.894	28.33	female
2943_AxiomGWHuSNP1_2010...	Pass	0.955	102.058	101.634	25.619	25.129	female
2944_AxiomGWHuSNP1_2010...	Pass	0.966	102.393	101.848	25.774	24.926	female
2947_AxiomGWHuSNP1_2010...	Pass	0.961	102.192	101.853	29.013	28.086	female
2948_AxiomGWHuSNP1_2010...	Pass	0.955	102.316	101.84	25.634	24.856	female
NAI1882_AxiomGWHuSNP1_...	Pass	0.915	101.553	101.186	28.796	27.976	female
NAI2057_a550094-4096756-0...	Pass	0.992	100.598	100.325	27.79	27.649	female
NAI2239_AxiomGWHuSNP1_...	Pass	0.88	100.634	100.195	29.26	28.962	female
NAI2249_AxiomGWHuSNP1_...	Pass	0.924	101.656	101.173	28.095	27.622	female

Find in Table | Row Count: 226 Selected: 0 | Show Filtered Only

Clearing an Individual Filter

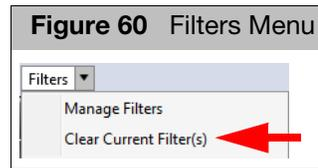
1. Right-click on the filtered column you want to clear.
The following window appears: (Figure 59)



2. Click **Clear Current Column Filter**.
The filter is removed.

Clearing All Current Filters

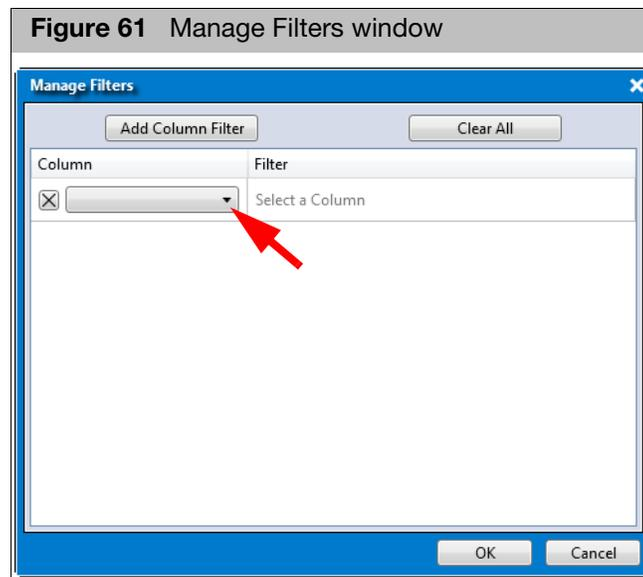
- Click the Filters drop-down, then select **Clear Current Filters**. (Figure 60)



Adding Filters (Method 2)

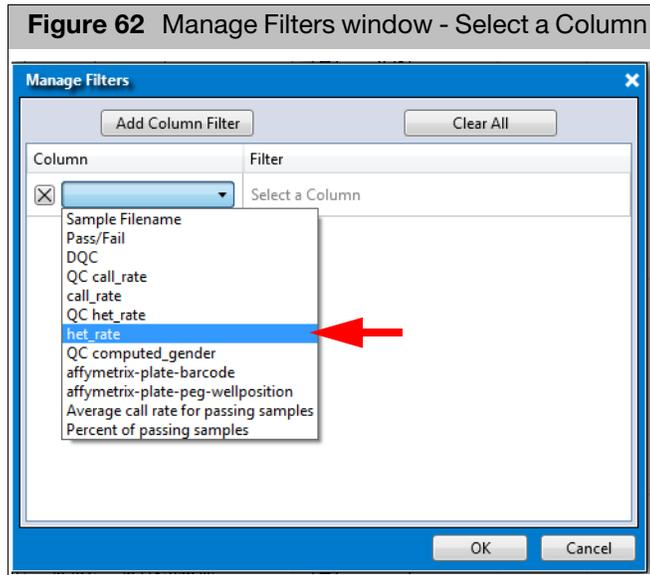
Use this method if you want to change more than one of your Sample Table column filters at the same time.

1. Click the **Filters** drop-down menu, then click **Manage Filters**.
The Manage Filters window appears. (Figure 61)

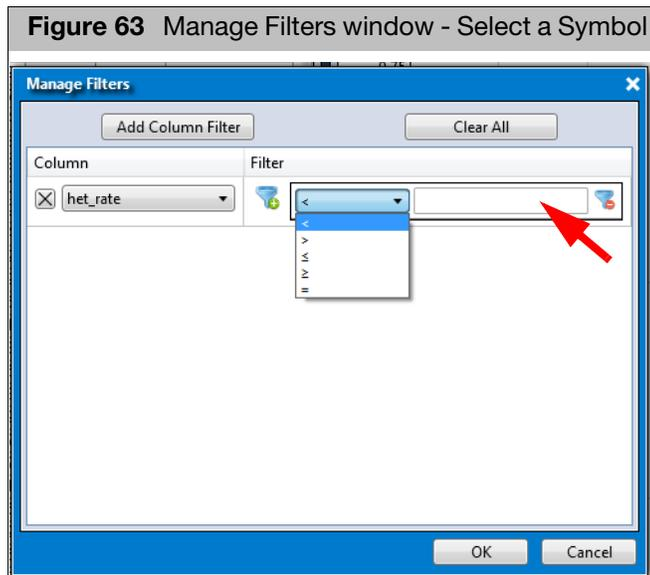


Note: If the column you want to filter contains text-based data, the Contains drop-down menu appears. If the column you want to filter contains numeric data, a symbol drop-down menu appears.

2. Click the **Column** drop-down, then click to select the Column name you want to create a filter for. (Figure 62)

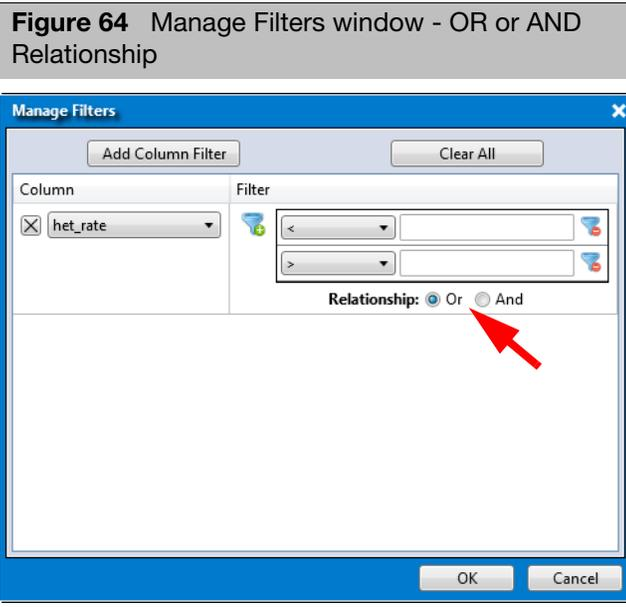


3. Click the **Symbol Value** drop-down menu to select the filtering symbol you want. (Figure 63)

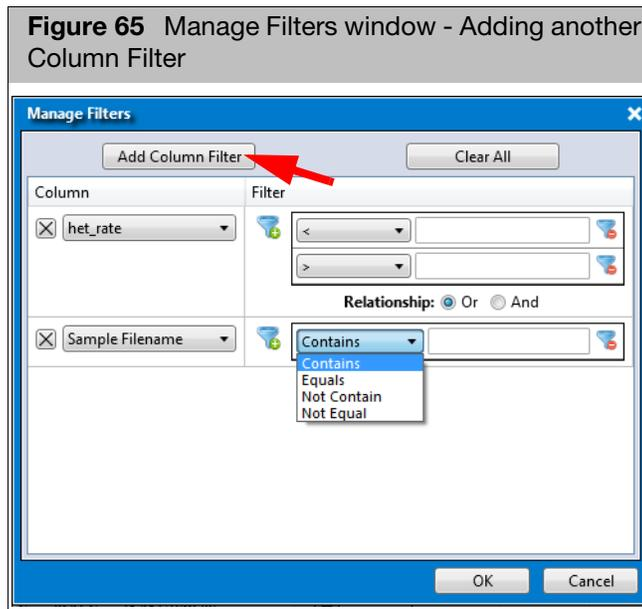


4. Click inside the text entry box to enter new value(s). (Figure 63)
5. OPTIONAL: If you want to add an additional filter to a column, click .

6. Click the **Or** or **And** radio button to choose **Or** or **AND** relationship logic. (Figure 64)



7. If needed, click **Add Column Filter**, then repeat the above steps. (Figure 65)

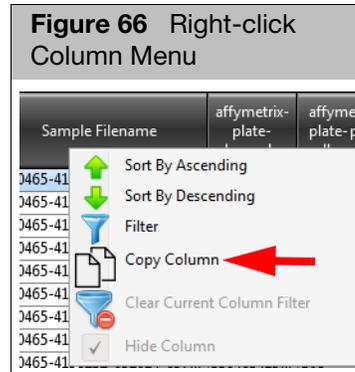


8. Click **OK**.
To remove a filter(s), click .
Click **Clear All** to remove ALL filters in the Manage Filters window.

Copying Column Data

To copy column data to your clipboard:

1. Click to select a column you want to copy to a clipboard, then right-click on it.
The following window appears: (Figure 66)



2. Click **Copy Column**.

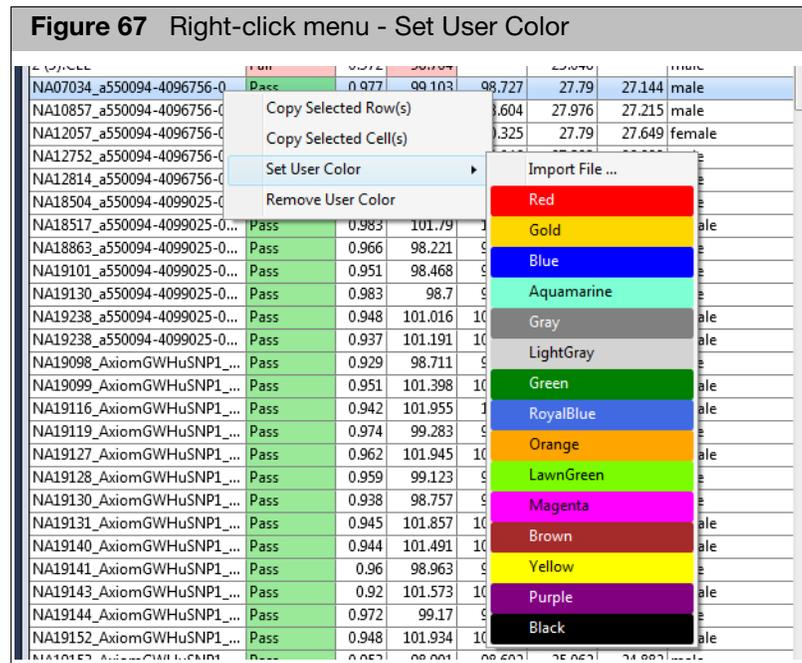
The column data is now ready for pasting (Ctrl v).

Setting User Colors

Use this feature to more easily identify different sets between the Sample Table and Cluster Graph.

Assigning a Color to a Sample

1. Right-click on the sample you want to assign a color to.
A menu appears. (Figure 67)

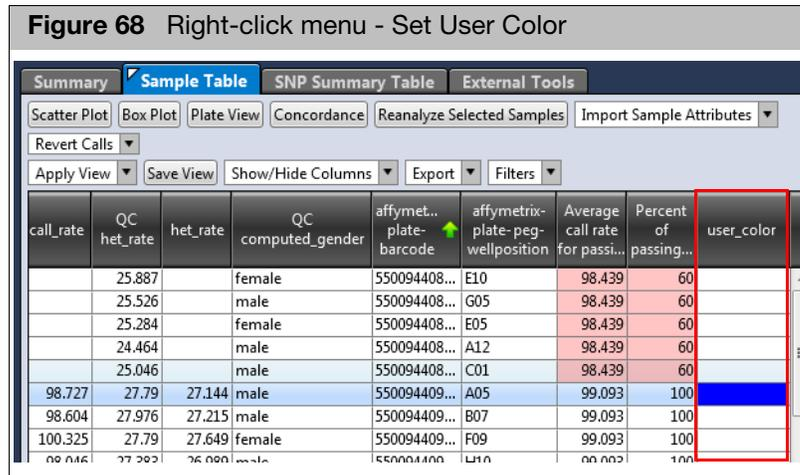


2. Mouse over **Set User Color**.

A color pallet appears.

- Click on the color you want.

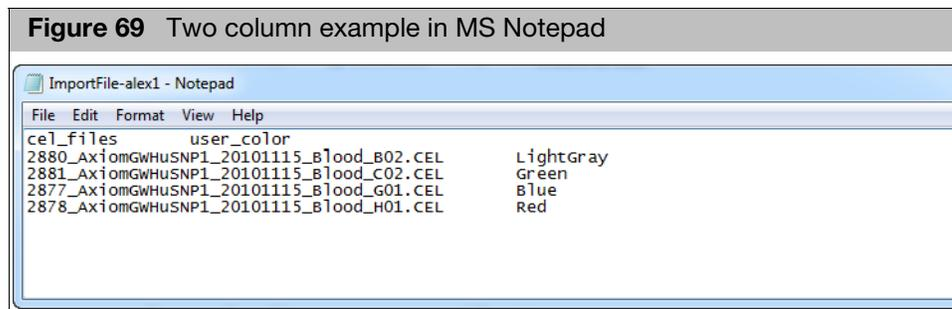
A **user_color** column is automatically added to your Sample Table and contains your sample's newly assigned color, as shown in [Figure 68](#).



Importing Assigned Colors

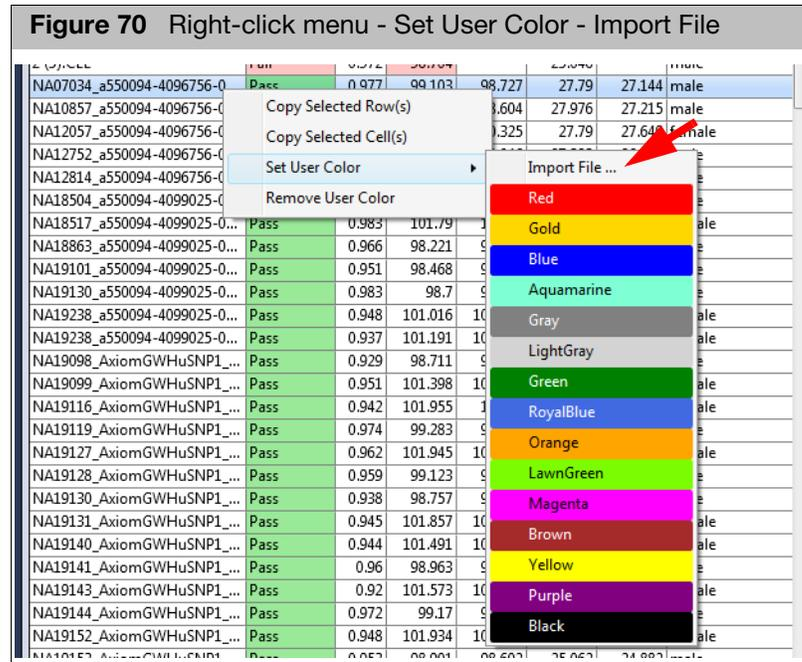
Use this feature if you want to assign colors to a large number of samples or if your Sample Table contains a vast amount of samples and you want to assign a color to only a few samples.

- Use MS Excel or MS Notepad (as you normally would) to create a two column table. ([Figure 69](#))



IMPORTANT! Your **user_color** entries must match the color pallet naming conventions shown in [Figure 70](#). Example: **RoyalBlue** not **Royal Blue**.

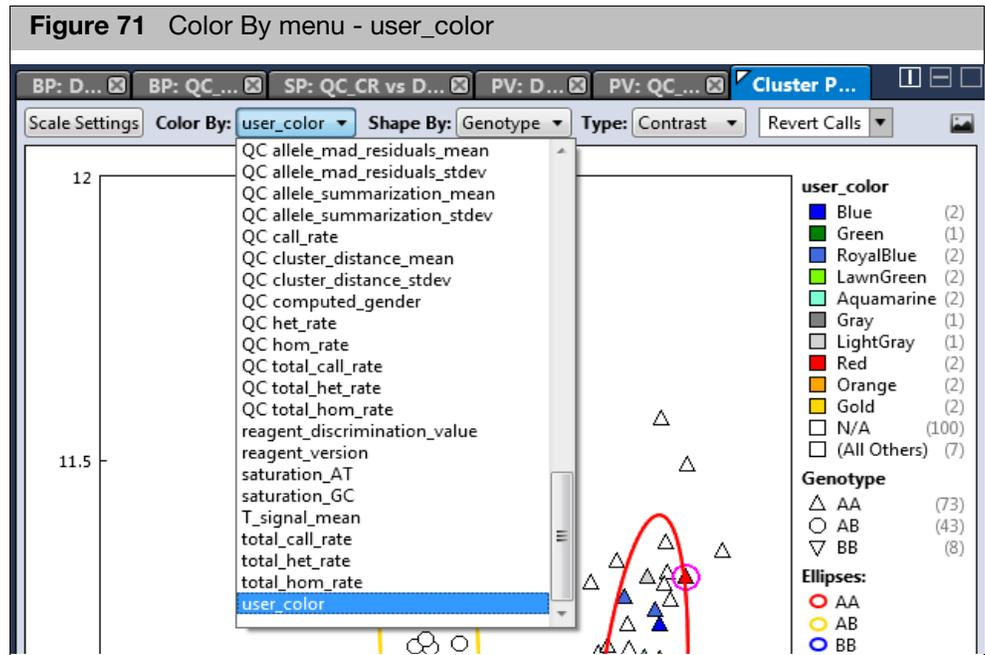
- Save your two column table as a tab-delimited text file to an easily accessible location.
- At the Sample Table, right-click on any sample.
A menu appears. ([Figure 70](#))



4. Mouse over **Set User Color**.
5. Click on **Import File...**
An **Import User Colors** Explorer window appears.
6. Locate your saved TXT file, click to highlight it, then click **Open**.
Your TXT file is now incorporated into the Sample Table.
7. Scroll the Sample Table right to see the added **user_color** column and assigned sample colors.

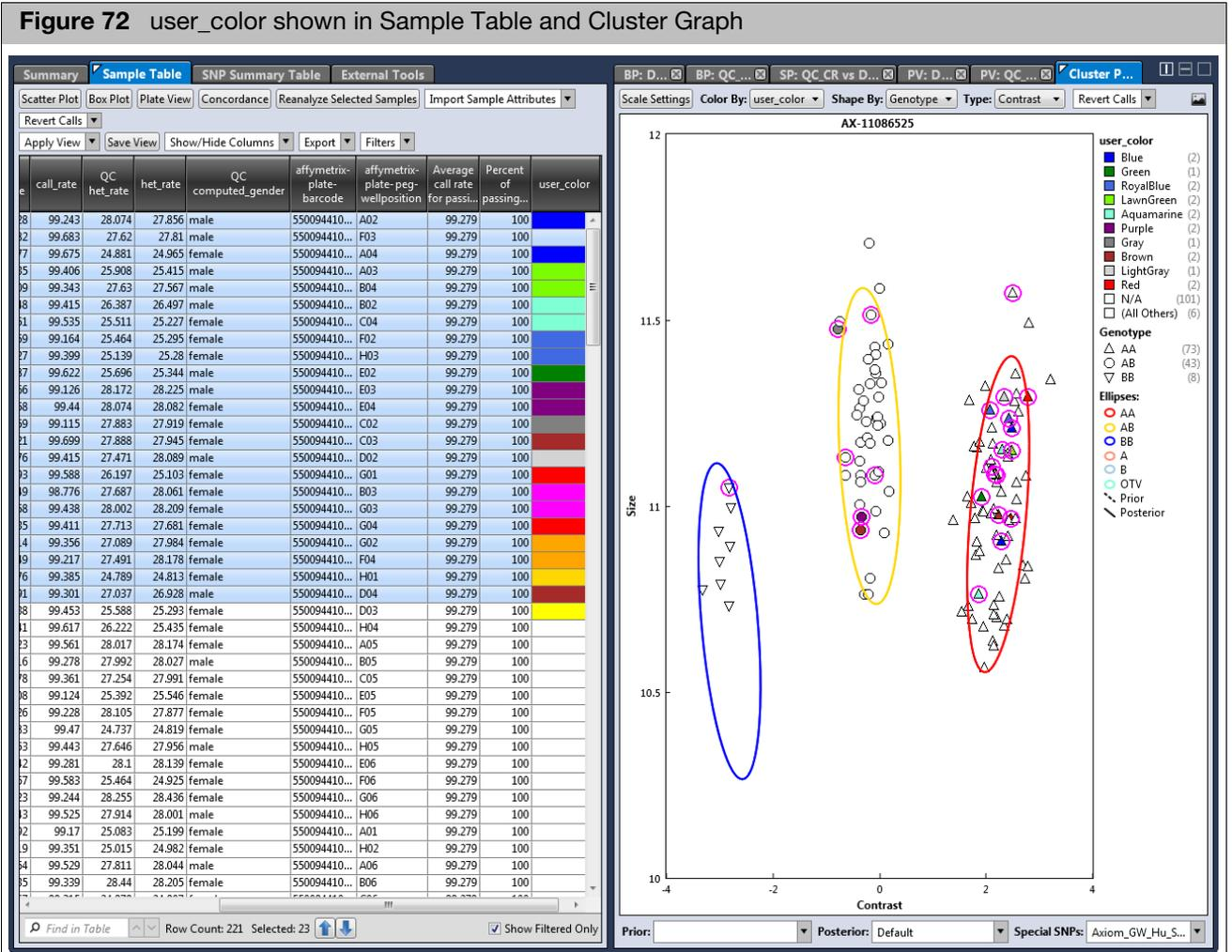
Viewing User Colors in the Cluster Graph

1. From the Cluster Graph, click the **Color By** drop-down menu. (Figure 71)



2. From the Sample Table, single-click on a color-coded sample file or Ctrl click, Shift click, or press Ctrl A (to select multiple color-coded sample files).

Each highlighted sample and its assigned color are now displayed in the Cluster Graph, as show in Figure 72.



Removing an Assigned User Color

1. From the Sample Table, right-click on the sample containing the color you want to remove.
A menu appears.
2. Click **Remove User Color**.
Your previously assigned sample color is now removed.

Searching Keywords

Note: The Find in Table tool can locate exact (case insensitive) matches. It also accepts wild-card (*) characters to aid in your search. Example: ABC*

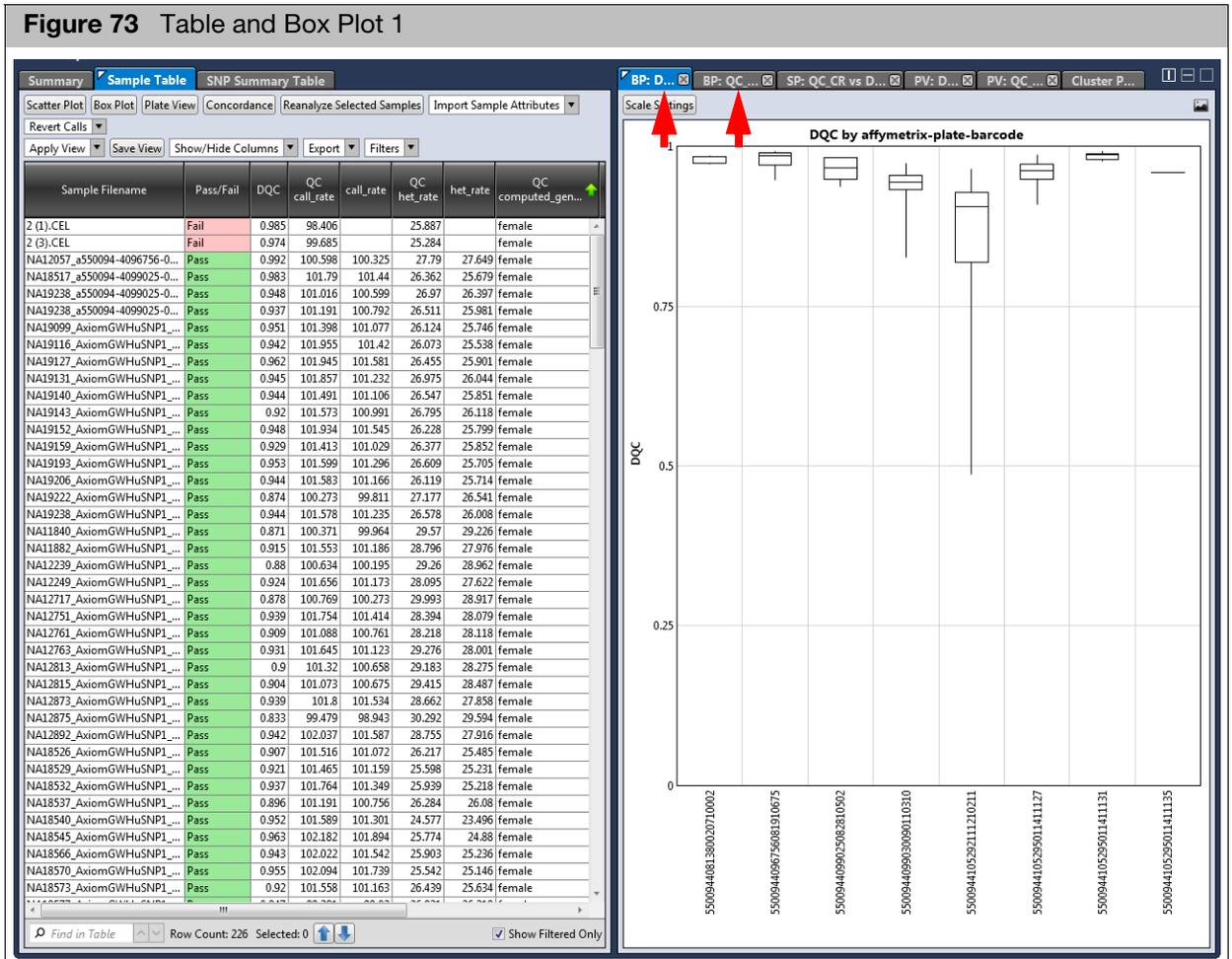
1. Click inside the  text field (bottom left corner of table).
2. Enter a keyword or number.
3. Click the **Up** or **Down** button.

When a match is found, the appropriate table entry is highlighted. If a graph is displayed, the appropriate graph point is also highlighted.

Box Plots

Note: By default, the Viewer generates two Box Plots. (Figure 73)

Viewing the Default Box Plots



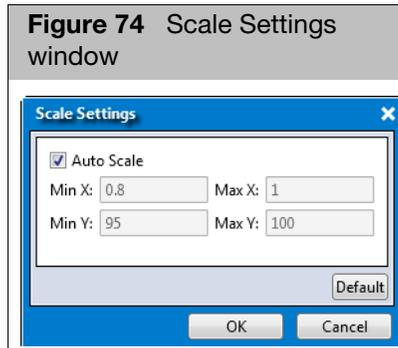
IMPORTANT! You cannot change a plot's axis values after it has been created. However, you can change its scale and coloring properties. See "Changing the Box Plot's Scale Setting Ranges" on page 62.

To change a Box Plot's axis properties, you must create a new Box Plot. See "Adding a New Box Plot" on page 62.

Changing the Box Plot's Scale Setting Ranges

1. Click **Scale Settings**.

The following window appears. (Figure 74)



By default, the window displays your current range values.

2. Uncheck the **Auto Scale** check box to enter different ranges in the provided text fields.
3. Click **OK**.

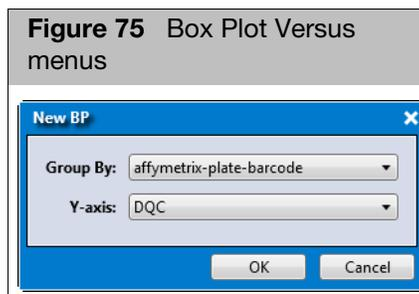
Your new settings are now reflected within the Box Plot. Modified Set Scale values are auto-saved.

If needed, click **Default** to return all values back to their factory settings.

Adding a New Box Plot

1. Click the **Box Plot** button.

The following window appears: (Figure 75)

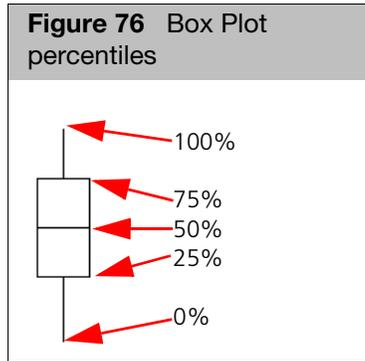


2. Click the **Group By** drop-down menu to select the X-axis for your new Box Plot. Your X-axis selection determines your new Box Plot's boxes and whiskers, based on the data group of values that are compiled.
3. Click the **Y-axis** drop-down menu to select the Y-axis you want. For Group By and Y-axis definitions, see "[Sample Table](#)" on page 154.
4. Click **OK**.

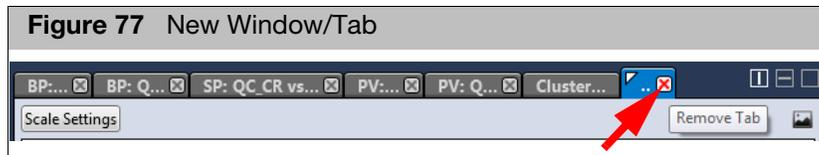
A new Box Plot window tab is created.

Reading Box Plot Percentiles

See [Figure 76](#).



At any time, click **X** to remove a window/tab, as shown in [Figure 77](#).



Saving the Current Box Plot View

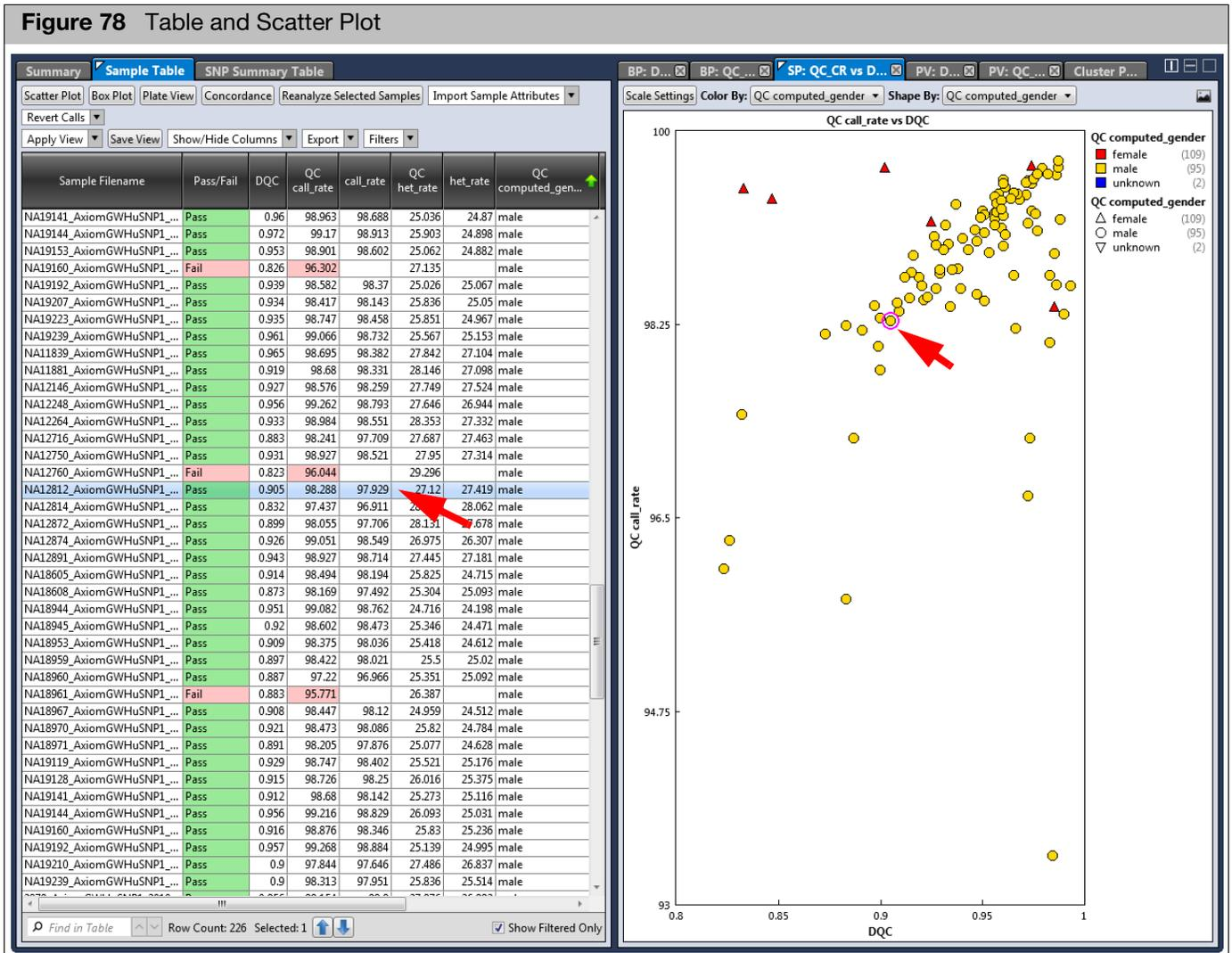
1. Click the **Save Image**  button.
An Explorer window appears.
2. Navigate to where you want to save the .PNG file, enter a filename, then click **Save**.

Scatter Plot

By default, the Viewer generates 1 Scatter Plot of QC call_rate vs. DQC. The data displayed in the plot are colored and shaped by QC computed_gender, as shown in Figure 78.

Viewing the Default Scatter Plot

1. Click to highlight a table entry to view its location within the Scatter Plot or click on a data point to highlight its corresponding entry in the Sample Table. (Figure 78)



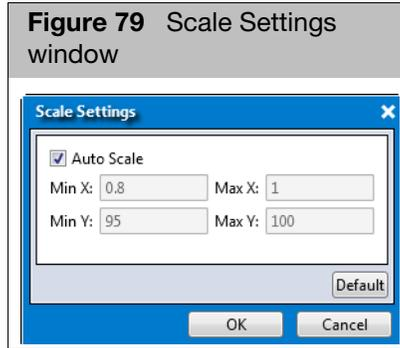
IMPORTANT! You cannot change the default Scatter Plot's pre-defined X and Y definitions, however you can change its Scale Settings and Color By and Shape By configuration.

To change a Scatter Plot's axis properties, you must create a new Scatter Plot. See "Adding a New Scatter Plot and Selecting its X and Y Properties" on page 65.

Changing the Scatter Plot's Setting Ranges

1. Click **Scale Settings**.

The following window appears. (Figure 79)



By default, the window displays your current range values.

2. Enter your new ranges in the appropriate text fields.
3. Click **OK**.

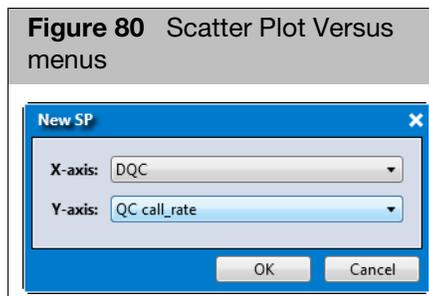
Your new settings are now reflected within the Scatter Plot. Modified Set scale values are auto-saved.

If needed, click **Default** to return all values back to their factory settings.

Adding a New Scatter Plot and Selecting its X and Y Properties

1. Click the **Scatter Plot** button.

The following window appears: (Figure 80)



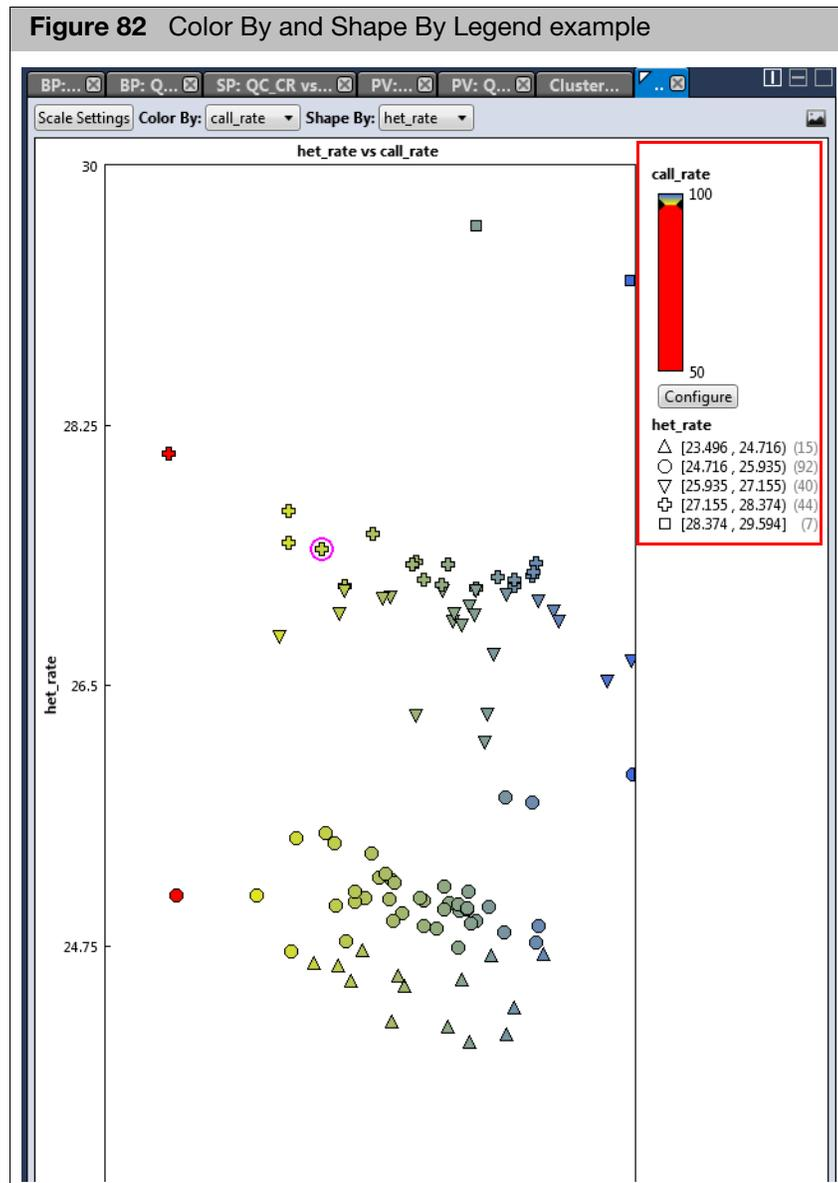
2. Use the drop-down menus to select your Plot's versus scenario (X and Y axis). For definitions, see "Sample Table" on page 154.
3. Click **OK**.

A new Scatter Plot window tab is created.

At any time, click **X** to remove a window/tab, as shown in Figure 81.



- Click the **Color By** and **Shape By** drop-down menus to select the combination view you want. See "Sample Table" on page 154 for Color By and Shape By definitions. Note: Your imported sample attributes are also available for use within the Color By and Shape By drop-down menus. For information on importing sample attributes, see "Importing Sample Attributes" on page 46. A legend appears within the plot. (Figure 82)



The graph can display up to 10 different colors and up to 10 different shapes. If the attributes selected for display have more than 10 categories, categories 1 through 9 are displayed normally, but categories 10 and higher get grouped together.

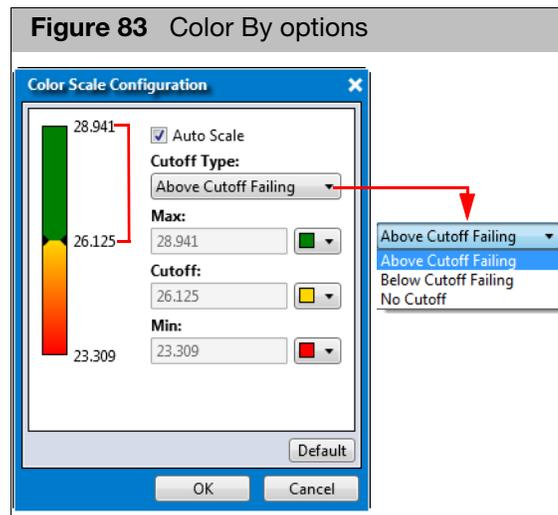
If your study has more than 10 values:

- If the value is text, the software takes the first nine values and assigns each a color or shape. The remaining values are put into a bin labeled **Other**. All values in the Other bin have the same color or shape.
- If the value is a date or number, the software divides the range of data into 10 equal bins and assigns a color or shape to each bin. If the data includes one or more outliers, it is possible to have one value in a particular bin and all other values in another bin.

Customizing Color By Settings

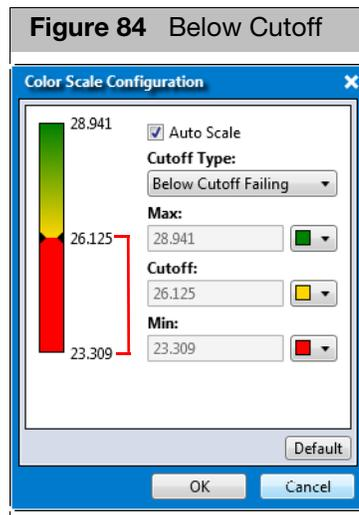
1. Click **Configure**.

The Color Scale Configuration window appears. (Figure 83)

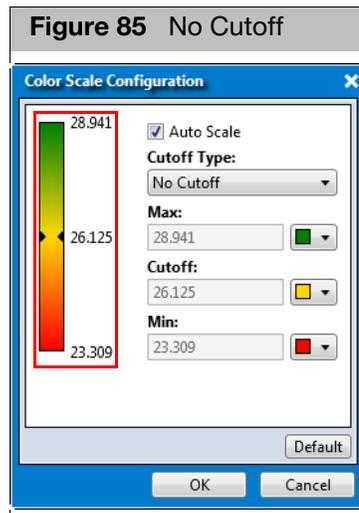


2. Use the provided text fields and color drop-down menus to customize your Color By selection.
 - **Auto Scale** check box (when checked) uses the actual minimum (lower bound) and maximum (upper bound) as your min/max scale. Uncheck the Auto Scale check box to enter your min and max number scales in the provided fields.
 - Click the **Cutoff Type** drop-down menu to select your cutoff preference.
 - **Above Cutoff Failing** - This presents a hard visual cutoff graph of all values that fail ABOVE the Cutoff value entered. The Above Cutoff data is represented by the color defined for Max. (Green in Figure 83)

- **Below Cutoff Failing** - This presents a hard visual cutoff graph of all values that fail BELOW the Cutoff value entered. The Below Cutoff data is represented by the color defined for Min. (Red in Figure 84)



- **No Cutoff** - This presents a smooth 3-point gradient of your defined Max, Min, and colors. (Figure 85)



3. Click **OK**.

Your Cutoff preference, entered values, and color selections are now displayed on the graph and saved for future use. If needed, click **Default** to revert all values back to their factory settings.

Saving the Current Scatter Plot View

1. Click the **Save Image**  button.
An Explorer window appears.
2. Navigate to where you want to save the .PNG file, enter a filename, then click **OK**.

Plate Views

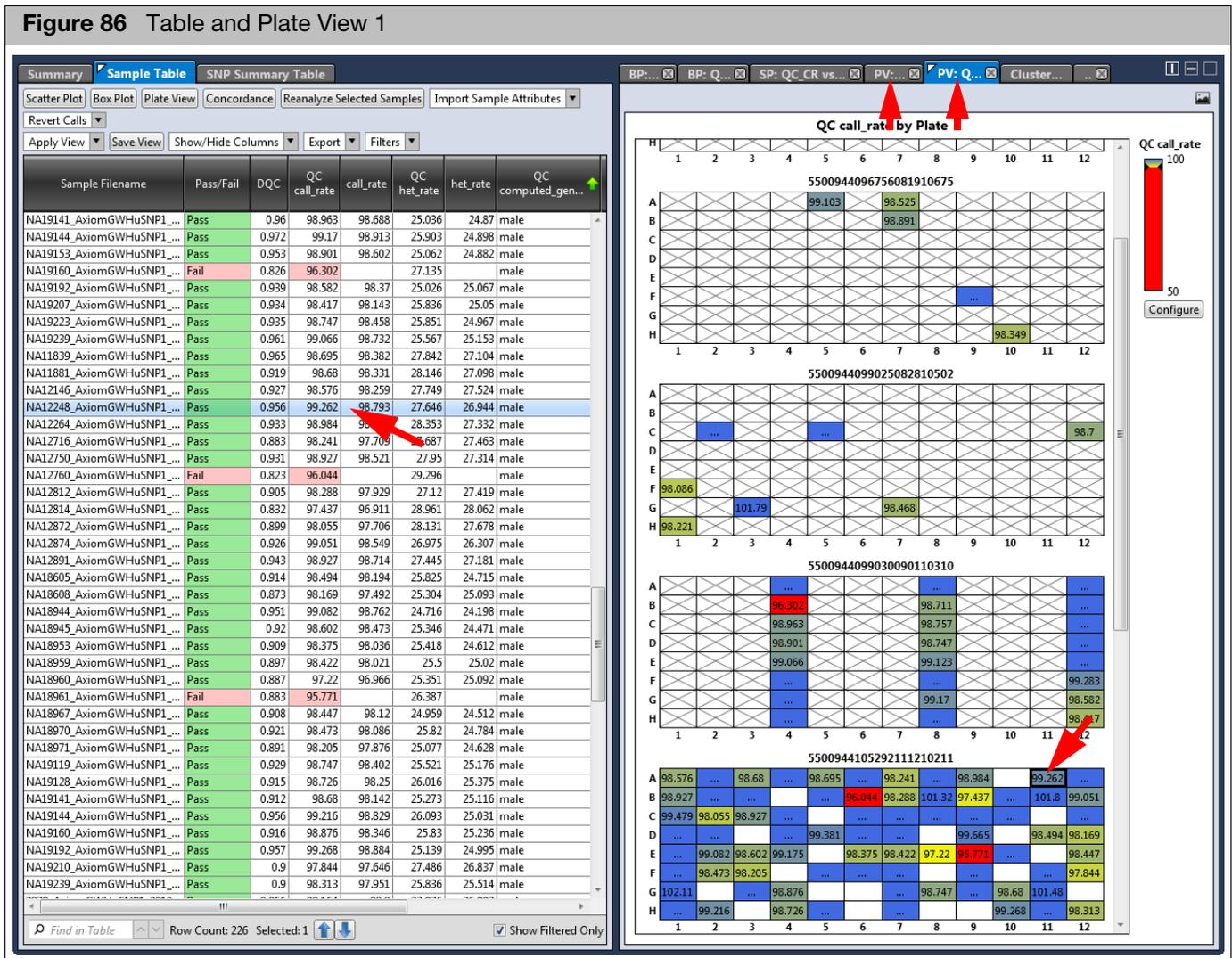
By default, the Viewer generates two Plate Views. (Figure 86) Each have a factory set metric. Plate View 1 is DQC by Plate and Plate View 2 is QC call_rate by Plate.

To display a different metric you must create a new Plate View. For more details, see "Adding a New Plate View Metric" on page 70.

The Plate Views display the currently selected (highlighted) metric from the Sample Table and are a graphic representation of the plate used. For example, 96 count plate layouts are shown in Figure 86.

Viewing the Default Plate Views

1. Click to highlight a table entry to view its location within the Plate View or click on a plate position to highlight its corresponding table entry, as shown in Figure 86.



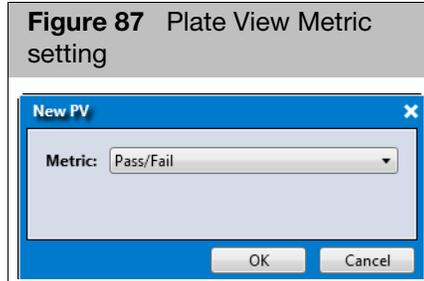
IMPORTANT! You cannot change a default Plate View, however you can change its Scale Settings, gradient, and coloring. See "Customizing your Plate View Settings" on page 71.

Adding a New Plate View Metric

The default Plate Views cannot be altered, therefore you must click the **Plate View** button to create a new Plate View to reflect your Metric change.

1. Click the **Plate View** button.

The following window appears: (Figure 87)



2. Use the drop-down menus to select your Plate View's Metric setting. See the tables in "Sample Table" on page 154 for Metric definitions.
3. Click **OK**.

The new Plate View window tab appears.

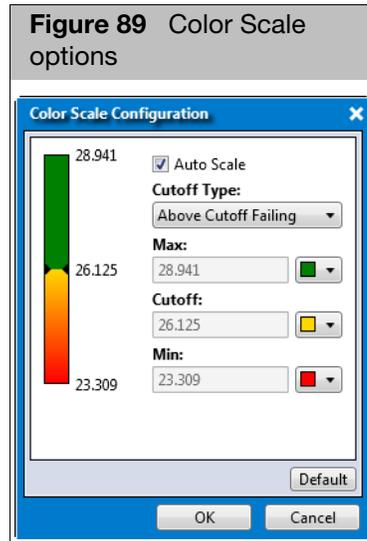
At any time, click **X** to remove a window/tab, as shown in Figure 88.



Customizing your Plate View Settings

1. Click **Configure** .

The Color Scale Configuration window appears. (Figure 89)



2. Use the provided text fields and color drop-down menus to customize your Color By selection.
 - **Auto Scale** check box (when checked) uses the actual minimum (lower bound) and maximum (upper bound) as your min/max scale. Uncheck the Auto Scale check box to enter your min and max number scales in the provided fields. Note: If the Auto Scale check box remains unchecked, you must enter new scale limits for each subsequent analysis.
 - Click the **Cutoff Type** drop-down menu to select the appropriate cutoff (based on the custom Cutoff value you entered).
3. Click **OK**.

Your new preferences are now displayed and saved for future use.

At any time, click the **Default** button to revert all the Color Scale Configuration window values back to their factory setting.

Saving the Current Plate View

1. Click the **Save Image**  button.

An Explorer window appears.

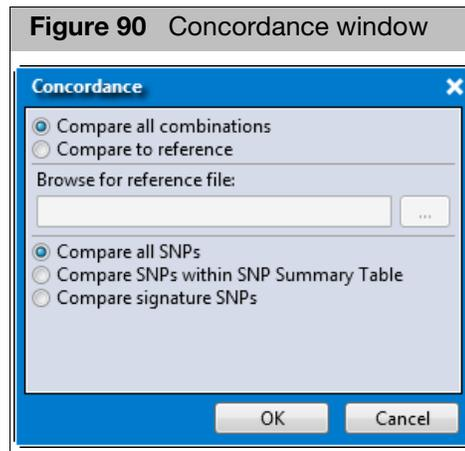
2. Navigate to where you want to save the .PNG file, enter a filename, then click **OK**.

Concordance Checks

- **Compare all combinations** enables you to compare the SNP calls for all samples. The concordance between all pairwise comparisons for the samples in the dataset/suitcase are reported.
- **Compare to reference** enables you to compare every sample to a single reference file.

Running a Concordance Check

1. Click the **Concordance** button.
The following window appears: (Figure 90)



Comparing All Combinations

1. Make sure the **Compare all combinations** radio button is selected.
2. By default, the **Compare all SNPs** button is selected. If needed, click the **Compare signature SNPs within the SNP Summary Table**, or **Compare signature SNPs** radio button.
3. Click **OK**.

After a few moments, the following Concordance window tab appears: (Figure 91)

X IMPORTANT! The amount of time to calculate concordance is proportional to the number of samples squared and the number of SNPs. It is highly recommended you use <1000 SNPs for an All versus All concordance check.

Figure 91 Default Concordance Table example

Index	Base	Reference	% Concordance
1	2877_Axio...	2878_Axio...	64.294
2	2877_Axio...	2879_Axio...	57.144
3	2877_Axio...	2880_Axio...	59.274
4	2877_Axio...	2881_Axio...	58.883
5	2877_Axio...	2882_Axio...	57.157
6	2877_Axio...	2883_Axio...	64.206
7	2877_Axio...	2884_Axio...	64.068
8	2877_Axio...	2885_Axio...	56.833
9	2877_Axio...	2887_Axio...	64.132
10	2877_Axio...	2888_Axio...	58.584
11	2877_Axio...	2889_Axio...	57.007
12	2877_Axio...	2890_Axio...	63.819
13	2877_Axio...	2891_Axio...	56.931
14	2877_Axio...	2892_Axio...	57.061
15	2877_Axio...	2893_Axio...	57.12
16	2877_Axio...	2894_Axio...	64.313
17	2877_Axio...	2895_Axio...	64.35
18	2877_Axio...	2896_Axio...	57.142
19	2877_Axio...	2897_Axio...	64.04
20	2877_Axio...	2898_Axio...	58.86
21	2877_Axio...	2899_Axio...	57.049
22	2877_Axio...	2900_Axio...	56.827
23	2877_Axio...	2901_Axio...	57.112
24	2877_Axio...	2902_Axio...	64.178
25	2877_Axio...	2903_Axio...	58.895
26	2877_Axio...	2904_Axio...	56.902
27	2877_Axio...	2905_Axio...	57.054
28	2877_Axio...	2907_Axio...	64.122
29	2877_Axio...	2908_Axio...	56.957
30	2877_Axio...	2909_Axio...	64.429
31	2877_Axio...	2910_Axio...	56.887
32	2877_Axio...	2915_Axio...	56.952
33	2877_Axio...	2916_Axio...	64.49
34	2877_Axio...	2917_Axio...	57.753
35	2877_Axio...	2918_Axio...	57.146
36	2877_Axio...	2943_Axio...	64.28
37	2877_Axio...	2944_Axio...	64.369
38	2877_Axio...	2946_Axio...	57.067
39	2877_Axio...	2947_Axio...	58.427
40	2877_Axio...	2948_Axio...	64.345

For definitions of the Concordance columns, see [Table 13](#) on page 158.

Comparing to Reference

1. Click the **Compare to reference** button.
The **Browse for the reference file** field is now activated.
2. Click the **Browse** button.
A Windows Explorer window appears.
3. Navigate to the appropriate reference file location, then click **Open**.
Your Reference file is displayed.

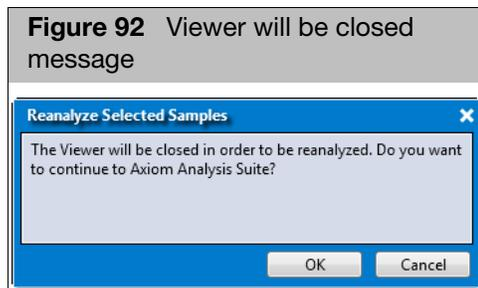
✗ IMPORTANT! The reference file you select can have more than two columns. However, only the first two columns are used during the *Compare to Reference* concordance check. Also, your reference genotypes must be reported using letter call codes (e.g. AA, AB, BB, NoCall).

4. By default, the **Compare all SNPs** button is selected. If needed, click the **Compare signature SNPs within the SNP Summary Table**, or **Compare signature SNPs** radio button.
5. Click **OK**.
After a few moments, a Concordance window tab appears.

Reanalyzing Samples

To reanalyze sample(s) displayed in the Sample Table:

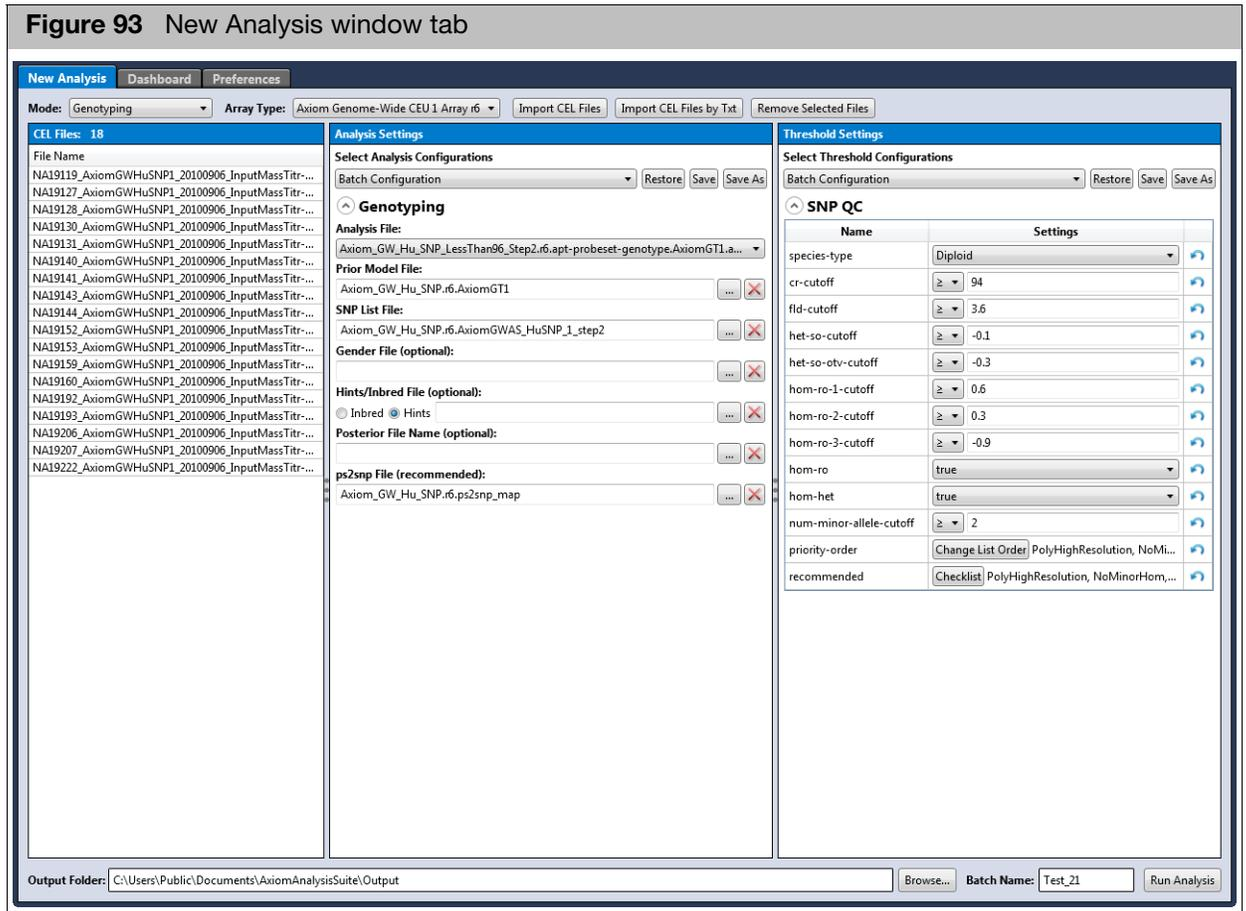
1. Single-click on a CEL file or Ctrl click, Shift click, or press Ctrl A (to select multiple files).
2. Click the **Reanalyze Selected Samples** button.
The following message appears. (Figure 92)



3. Click **OK**.

The Viewer closes. Your selected samples are displayed and ready for re-analysis in a New Analysis window tab. (Figure 93)

Figure 93 New Analysis window tab



- See Chapter 2, "Performing an Analysis" on page 19 for instructions on setting up an analysis.

4

The Viewer: SNP Summary Table and Cluster Plot

SNP Summary Table

Figure 94 is an example of a standard SNP Summary Table

Figure 94 SNP Summary Table window

probeset_id	ConversionType	CR	MinorAlleleFrequency	H.W.p-Value	FLD	HomFLD	HetSO	HomRO	Nclus	n_AA	n_AB	n_BB	n_NC	hemizygous	HomHet	gender_metrics	Call Modified
AFFX-KIT-000001	PolyHighResol...	100	0.464	0.306	20.691	44.561	0.325	2.754	3	39	76	51	0	0	0	all	False
AFFX-KIT-000002	PolyHighResol...	99.398	0.33	0.482	8.127	18.035	0.191	0.846	3	76	69	20	1	0	0	all	False
AFFX-KIT-000003	PolyHighResol...	100	0.22	0.007	6.492	13.807	0.146	0.738	3	107	45	14	0	0	0	all	False
AFFX-KIT-000004	PolyHighResol...	100	0.319	0.145	15.227	33.186	0.318	2.124	3	81	64	21	0	0	0	all	False
AFFX-KIT-000005	PolyHighResol...	100	0.081	1	5.139	11.779	0.094	1.069	3	140	25	1	0	0	0	all	False
AFFX-KIT-000008	PolyHighResol...	100	0.322	0.532	5.412	11.248	0.133	1.127	3	19	69	78	0	0	0	all	False
AFFX-KIT-000009	NoMinorHom	100	0.075	0.602	21.081		0.247	4.47	2	0	25	141	0	0	1	all	False
AFFX-KIT-000012	PolyHighResol...	100	0.368	0.388	20.114	42.061	0.238	2.784	3	69	72	25	0	0	0	all	False
AFFX-KIT-000013	PolyHighResol...	99.398	0.136	0.19	5.134	10.814	0.169	0.813	3	125	35	5	1	0	0	all	False
AFFX-KIT-000014	PolyHighResol...	100	0.099	0.663	18.947	38.956	0.567	2.803	3	2	29	135	0	0	0	all	False
AFFX-KIT-000015	PolyHighResol...	100	0.181	1	30.972	67.77	0.439	4.658	3	5	50	111	0	0	0	all	False
AFFX-KIT-000016	PolyHighResol...	100	0.135	1	23.086	48.374	0.579	2.409	3	3	39	124	0	0	0	all	False
AFFX-KIT-000017	PolyHighResol...	100	0.078	1	20.903	44.214	0.42	3.888	3	141	24	1	0	0	0	all	False
AFFX-KIT-000018	PolyHighResol...	99.398	0.23	1	16.629	34.399	0.254	2.451	3	9	58	98	1	0	0	all	False
AFFX-KIT-000019	PolyHighResol...	99.398	0.373	0.043	6.281	15.899	0.235	1.536	3	29	65	71	1	0	0	all	False
AFFX-KIT-000021	PolyHighResol...	99.398	0.179	0.79	8.169	17.256	0.331	1.384	3	112	47	6	1	0	0	all	False
AFFX-KIT-000022	PolyHighResol...	100	0.434	0.381	18.57	38.557	0.397	2.523	3	56	76	34	0	0	0	all	False
AFFX-KIT-000023	PolyHighResol...	99.398	0.309	0.238	6.384	14.553	0.193	1.015	3	82	64	19	1	0	0	all	False
AFFX-KIT-000025	PolyHighResol...	100	0.494	0.439	17.222	36.978	0.348	2.333	3	43	78	45	0	0	0	all	False
AFFX-KIT-000026	PolyHighResol...	99.398	0.436	0.147	13.408	33.126	0.509	1.785	3	36	72	57	1	0	0	all	False
AFFX-KIT-000027	PolyHighResol...	100	0.41	0	23.078	47.688	0.238	2.894	3	41	54	71	0	0	0	all	False
AFFX-KIT-000029	PolyHighResol...	99.398	0.397	0	11.659	24.178	0.135	1.097	3	75	49	41	1	0	0	all	False
AFFX-KIT-000031	PolyHighResol...	100	0.238	0.671	15.932	33.638	0.379	2.353	3	95	63	8	0	0	0	all	False
AFFX-KIT-000032	PolyHighResol...	100	0.377	0.071	20.768	43.819	0.607	2.467	3	29	67	70	0	0	0	all	False
AFFX-KIT-000033	PolyHighResol...	99.398	0.258	0	13.454	27.084	0.505	2.629	3	102	41	22	1	0	0	all	False
AFFX-KIT-000049	PolyHighResol...	100	0.386	0.583	12.018	25.089	0.511	2.074	3	61	82	23	0	0	0	all	False
AFFX-KIT-000050	PolyHighResol...	100	0.473	0.227	6.032	12.798	0.147	0.776	3	50	75	41	0	0	0	all	False
AFFX-KIT-000121	PolyHighResol...	100	0.283	0.15	13.623	28.574	0.309	2.419	3	10	31	49	0	0	0	female	False
AFFX-KIT-000198	PolyHighResol...	100	0.356	0.228	17.117	36.866	0.365	2.704	3	40	36	14	0	0	0	female	False
AFFX-KIT-000209	PolyHighResol...	100	0.333	0.018	4.766	10.896	0.129	0.697	3	15	30	45	0	0	0	female	False
AFFX-KIT-000226	PolyHighResol...	98.889	0.371	0.002	12.302	27.039	0.556	2.075	3	42	28	19	1	0	0	female	False
AFFX-KIT-000408	PolyHighResol...	100	0.217	1	19.59	41.485	0.12	3.567	3	4	31	55	0	0	0	female	False
AX-162003418	MonoHighRes...	100								1						NA	False
AX-162005580	PolyHighResol...	100								2						NA	False
AX-162007660	CallRateBelow...	77.711								3						NA	False
AX-162011899	PolyHighResol...	100								3						NA	False
AX-162020167	Other	100								1						NA	False

NA Row Count: 37 Selected: 0 Show Filtered Only

Multi-allele probesets are probesets that can report calls from more than two alleles. Multi allele probesets can be identified by selecting the annotation field **Allele_Count**. Additional columns can be shown that report metrics for multi-allele probesets.

Multi-allele probesets are reported differently than the more common biallele probesets.

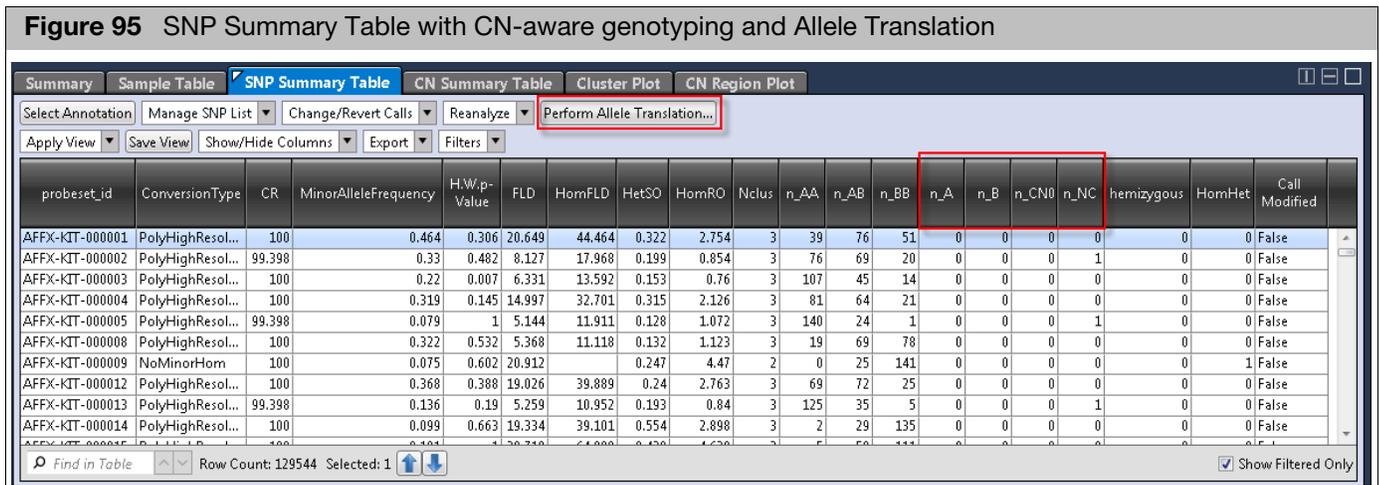
- Multi-allele probesets will not report NoMinorHom or OTV ConversionTypes.
- Multi-allele probesets do not report values for many of the default SNP Summary Table columns. Click **Show/Hide Columns** to select additional Columns with multi-allele probeset metrics.

Note: If an analysis batch does not include samples needed to evaluate a SNP's performance, then no metrics (including ConversionType) will be reported. For example, if an analysis batch does not contain male samples, no metrics are reported for Y chromosome SNPs and none of these SNPs are added to the Recommended SNP list.

Figure 95 represents a SNP Summary Table with CN-aware genotyping and Allele Translation.

Note: The **Perform Allele Translation...** button only appears if supported arrays are available. See Chapter 5, "Allele Translation" on page 97 for more information.

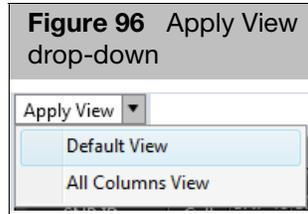
A SNP Summary Table with CN-aware genotyping also displays a count of calls for hemizygous genotypes (n_A, n_B) and a count of ZeroCN calls (n_CN0), as shown in Figure 95. These additional columns appear for arrays that support copy number-aware genotyping. For more information, see Appendix A, "Predefined Region CN Analysis Variations" on page 122.



Using the SNP Summary Table

Setting your SNP Summary Table View

1. Click the **Apply View** drop-down. (Figure 96)

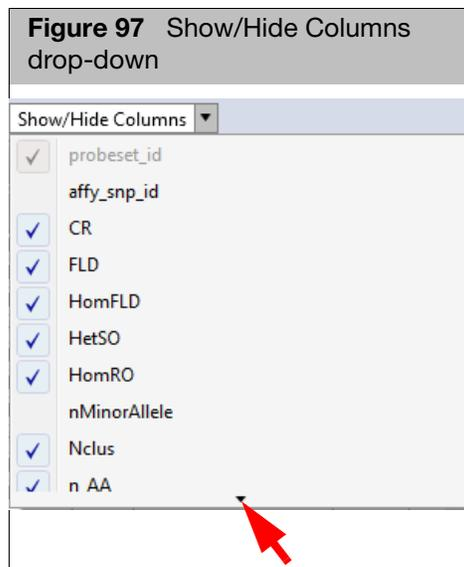


- The **Default View** is the initial table view and includes a preset number of columns.
- The **All Columns View** displays the maximum available columns.

Adding and Removing Table Columns

1. Click the **Show/Hide Columns** drop-down. (Figure 97)

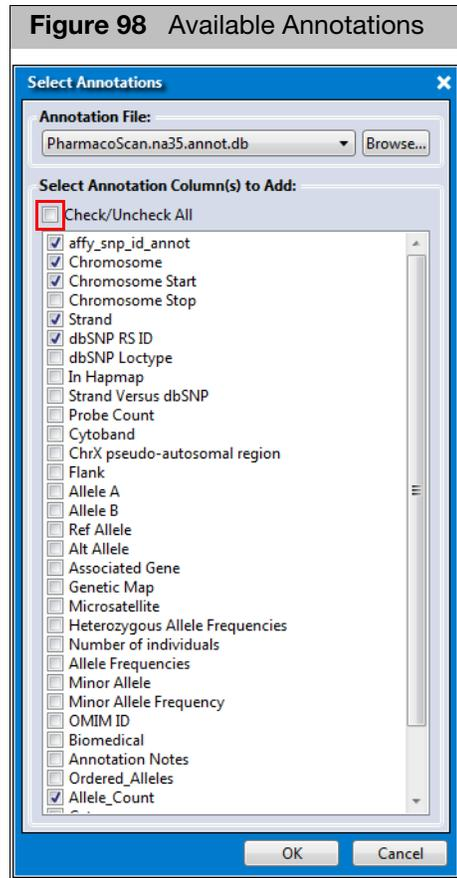
A list of available columns appear.



2. Click the check box next to the column(s) you want to add (Show) on the table. Click to uncheck a column you want to remove (Hide) from the table. Mouse over the menu's down arrow to reveal more available column choices. See [Table 15 on page 162](#) for column definitions.

Selecting Annotations

1. Click the **Select Annotation** button.
The following window appears. (Figure 98)



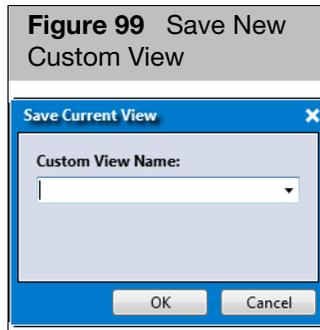
2. Select the appropriate Annotation File from the drop-down menu list.
3. Click the check box next to the Annotation Column(s) you want to add to the table or click the **Check/Uncheck All** check box (Figure 98) to add or remove ALL available annotations. See "[Annotations and Columns](#)" on page 159 for each Annotation's definition.
4. Click **OK**.

Your selected annotation columns are now added to the right side of the SNP Summary Table.

Saving your Table Column View

1. After you have your preferred SNP Summary Table columns set, click **Save View**.

The following window appears: (Figure 99)



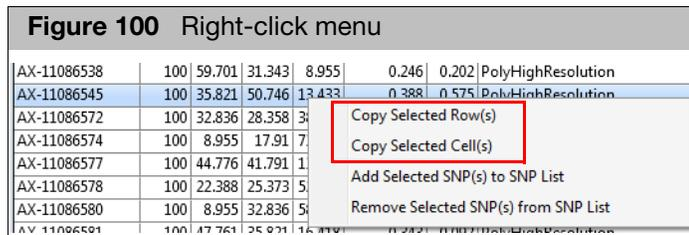
2. Enter a name, then click **OK**.

Your custom table view is now saved and stored inside the Apply View menu for future use.

Copying Selected Row(s)

1. Right-click on a row you want to copy. Optional: Single-click on a row or Ctrl click, Shift click, or press Ctrl A (to select multiple rows).

A menu appears. (Figure 100)



2. Click **Copy Selected Row(s)**.

The row data is now ready for pasting (Ctrl v).

Copying Selected Cell(s)

1. Right-click on a cell you want to copy. Optional: Single-click on a row or Ctrl click, Shift click, or press Ctrl A (to select multiple cells).

A menu appears. (Figure 100)

2. Click **Copy Selected Cell(s)**.

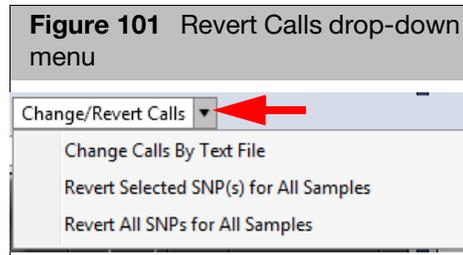
The cell data is now ready for pasting (Ctrl v).

Changing or Reverting Genotype Calls

Reverting Calls changes them back to what they were originally called by the algorithm. No other history is saved, only its current and original values.

1. Click the **Change/Revert Calls** drop-down.

The following menu appears: (Figure 101)



Changing Genotype Calls by Text File

1. Click **Change Calls By Text File**.

An Explorer window appears.

2. Navigate to the text file's location.

✗ IMPORTANT! Your Change Call by Text file must start with the header *probeset_id* and use numeric or alphabetic codes for the genotype, as shown in Figure 102. It also cannot contain any annotation columns.

Figure 102 Change Call by Text file example shown in Excel (as a tab-delimited text file)

	A	B	C	D	E	F	G	H	I	J	K	L	M	N	O	P
1	probeset_id	2877_Axic	2878_Axic	2879_Axic	2880_Axic	2881_Axic	2882_Axic	2883_Axic	2884_Axic	2885_Axic	2887_Axic	2888_Axic	2889_Axic	2890_Axic	2891_Axic	2892_A
2	AX-11699985	AA	AA													
3	AX-11700086	AA	AA													
4	AX-11700405	BB	BB	AB	BB	AB	AA	BB	BB	BB	AB	BB	AA	BB	AA	AA
5	AX-11700418	AB	AB	BB	BB	AB	AB	AB	BB	BB	BB	AB	BB	AB	BB	BB
6	AX-11700675	BB	BB													
7	AX-11700679	AB	BB	BB												
8	AX-11700939	AA	AA													

3. Click **Open**.

Your Genotype Calls are now changed and reflected in the Cluster Plot.

Reverting Selected SNPs for all Samples

1. Single-click on a SNP file or Ctrl click, Shift click, or press Ctrl A (to select multiple files) within the SNP Sample Table.
2. Click **Revert Selected SNP(s) for All Samples**.

Reverting all SNPs for all Samples

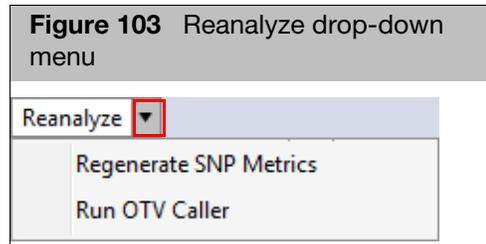
Use this feature to perform a master SNPs reset.

1. Click **Revert All SNPs for All Samples**.

✗ IMPORTANT! Once *Revert Calls* is performed, the selected calls will be reverted to original calls. This cannot be undone.

Reanalyzing your SNP Summary Table Data

The **Reanalyze** drop-down menu (Figure 103) offers optional steps for post-processing SNP data. These functions utilize the output files from the previous genotyping and classification steps as input.

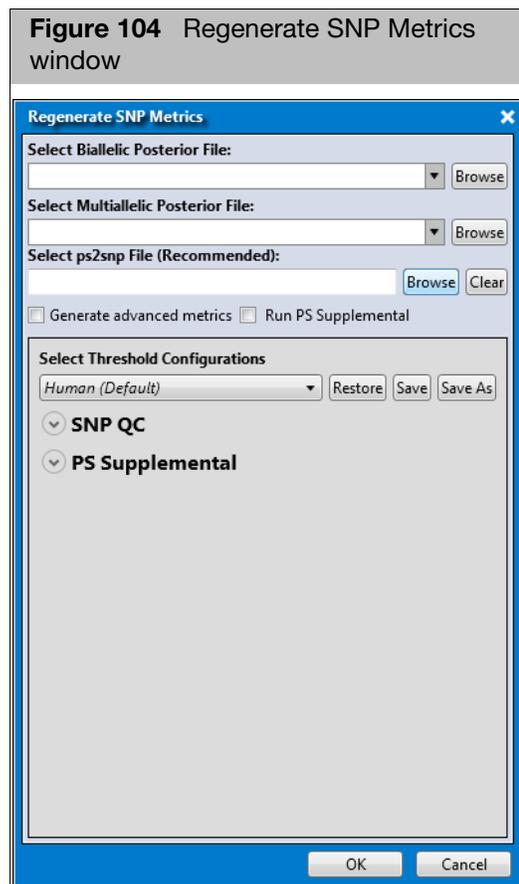


Regenerate SNP Metrics

The **Regenerate SNP Metrics** operation allows you to recalculate SNP Metrics and ConversionType classifications, which you may want to do if you have edited any genotype calls, or if you want to change SNP QC thresholds, or if you want to generate additional metrics and classifications.

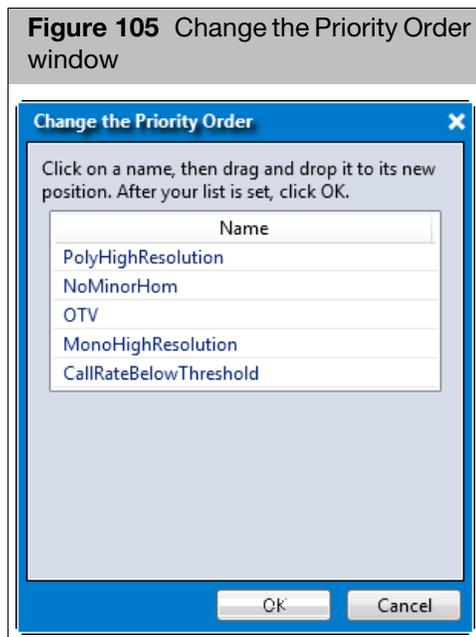
1. Click **Regenerate SNP Metrics**.

The following window appears: (Figure 104)



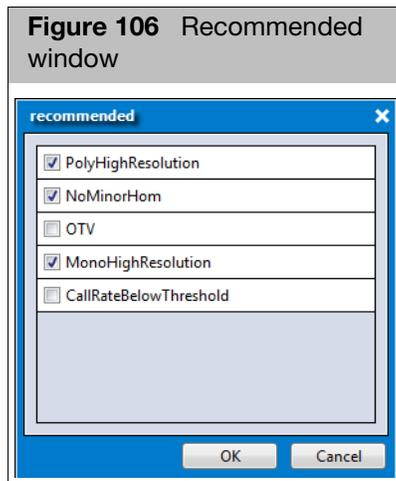
1. Use the drop-down to select the appropriate Posterior File or click its **Browse** button.
An Explorer window appears.
2. Navigate to Posterior File you want to use, then click **Open**.
3. Choose an appropriate ps2snp File (recommended), as described in "Using the Analysis Settings Fields" on page 24.
4. (Optional) Click the **Generate advanced metrics** check box to calculate additional SNP-specific posterior metrics.
5. (Optional) Click the **Run PS Supplemental** check box to generate additional metrics and probeset ConversionTypes. Note: PS Supplemental performs further classification that may be needed for polyploid organisms, complex genomes, or inbred populations.
6. Select the Threshold Configuration you want to use, as described in "Customizing Thresholds" on page 28.
7. Click to expand and use the **SNP QC** drop-down menu selections and text fields to setup the regeneration of your SNP Metric.
8. To change the priority-order of the various assigned ConversionTypes, go to the SNP QC setting's priority-order column and click the **Change List Order** button.

The following window appears: (Figure 105)



9. Click and hold onto the selection you want to move, then drag and drop it into its new position. After you get the order of priority you want, click **OK**.
10. To change the recommended options, click **Checklist**.

The following window appears: (Figure 106)



11. Click to check/uncheck the available recommended options. Click a field's  button to return its value back to its default setting.

Note: If you have checked **Run PS Supplemental** check box, then the associated threshold configurations are relevant. Edit them as needed.

12. Click **OK**.

After the process is complete, your SNP Summary Table is updated.

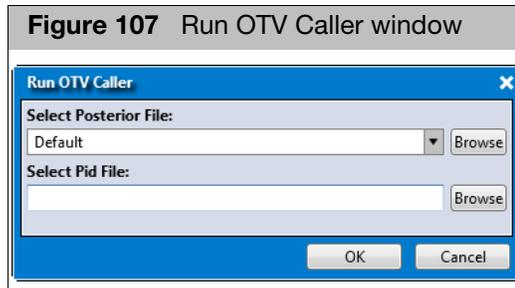
Running OTV Caller

The OTV Caller is intended for SNPs that have been classified as likely having off-target variants (OTV), or for SNPs with unusually large Y-dimension variance (as identified by PS Supplemental option of the Regenerate SNP Metrics).

OTV Caller function performs post-processing analysis to identify miscalled clustering and identify which samples should be in the OTV cluster and which samples should remain in the AA, AB, or BB clusters. Samples in the OTV cluster are re-labeled as OTV.

1. Click **Run OTV Caller**.

The following window appears: (Figure 107)



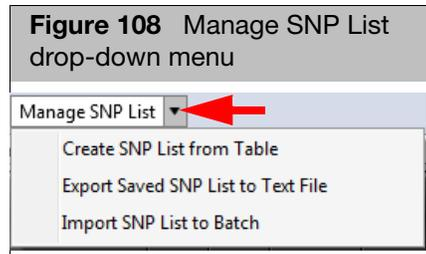
1. Use the drop-down to select the appropriate Posterior File or click its **Browse** button.
An Explorer window appears.
2. Navigate to Posterior File you want to use, then click **Open**.
3. Click the Select Pid File's **Browse** button.
An Explorer window appears.
4. Navigate to Pid File you want to use, then click **Open**.
5. Click **OK**.

Managing your SNP List

Axiom Analysis Suite enables lists of SNPs to be saved within the application.

1. Click the Manage SNP List drop-down menu (Figure 108), then select the option you want.

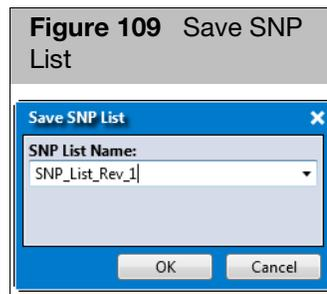
Note: The Recommended SNP List is auto-generated and updated whenever SNP metrics are calculated. Avoid using this reserved (default) SNP List name for your custom lists.



Saving your Current SNP List

1. To save all SNPs currently displayed in the SNP Summary Table, click **Create SNP List from Table**.

The following window appears: (Figure 109)



2. Enter a name, then click **OK**.

Exporting your SNP List

Before exporting a SNP List you must first create one. If no SNP Lists are detected, a message box appears. Click **OK** to acknowledge the message, then go to "Saving your Current SNP List" to create a SNP List.

1. Click **Export Saved SNP List to Text File**.

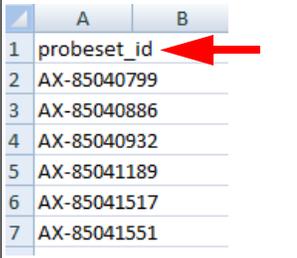
An Explorer window appears.

2. Navigate to an export location, enter a name, then click **Save**.

Importing a SNP List

1. Click **Import SNP List to Batch**.
An Explorer window appears.
2. Navigate to your SNP List location containing your tab-delimited text file.
Your first row/column header must be labeled **probeset_id**, as shown in [Figure 110](#), otherwise an error message appears.

Figure 110 Required probeset_id column header



	A	B
1	probeset_id	
2	AX-85040799	
3	AX-85040886	
4	AX-85040932	
5	AX-85041189	
6	AX-85041517	
7	AX-85041551	

3. Click **Open**.
Your imported SNP List now appears in the SNP Summary Table.

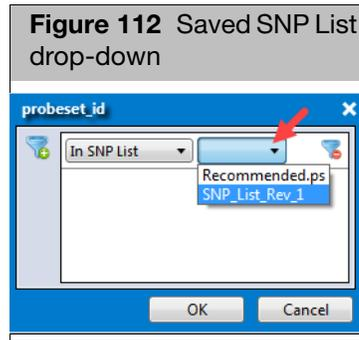
Using your Saved SNP List

Displaying SNPs in a SNP list

1. Click to select the **probeset_id** column, then right-click on it.
2. Click **Filter**.
The Filter window appears. ([Figure 111](#))



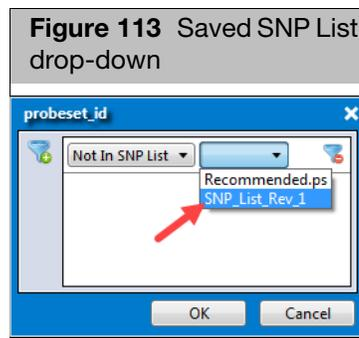
3. Add your previously saved SNP List by selecting it from the drop-down list. (Figure 112)



4. Click **OK**.
Only the SNPs in your SNP List are displayed in the SNP Summary Table.

Displaying SNPs that are not in your SNP List

1. Click to select the **probeset_id** column, then right-click on it.
2. Click **Filter**.
3. Click the **In SNP List** drop-down, then click to select **Not in SNP List**. (Figure 113)
4. Click the Saved SNP List drop-down, select your saved SNP List, then click **OK**.



Your SNPs from the SNP List are no longer displayed in the SNP Summary Table.

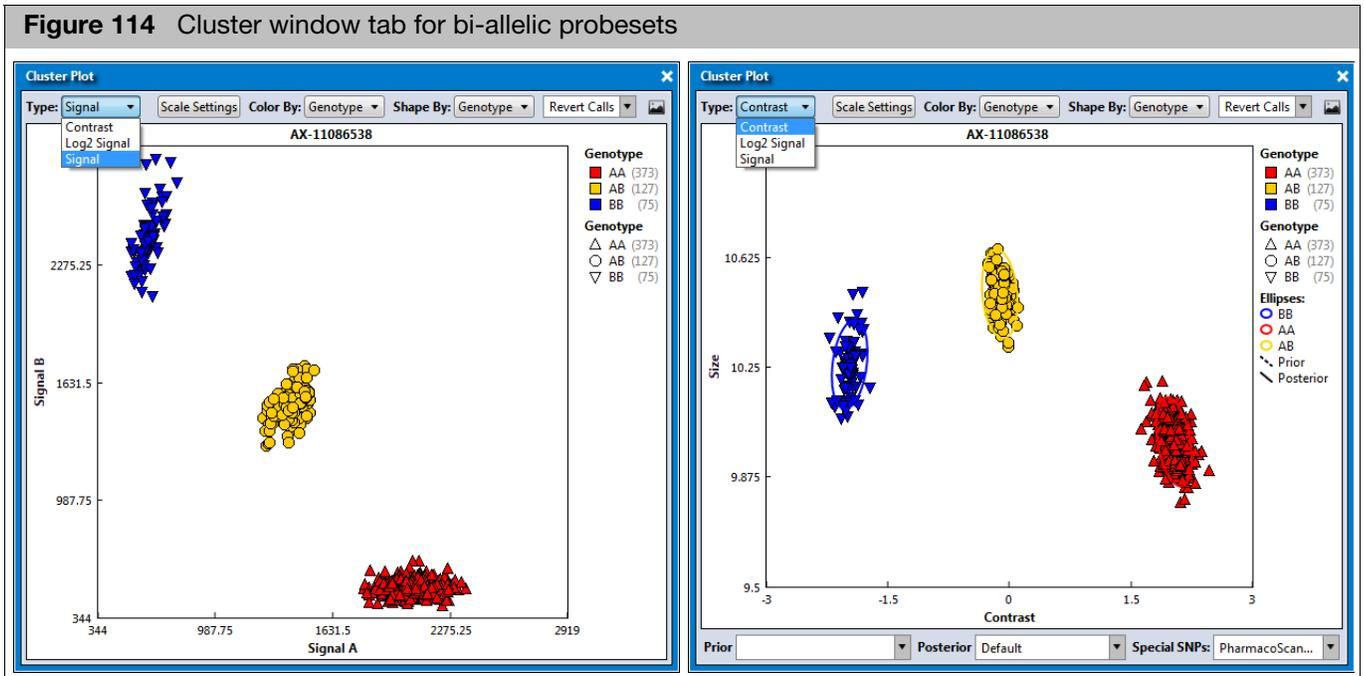
Cluster Plot

The Cluster Plot displays the SNP calls for selected samples as a set of points in the clustering space used for making the calls. A visual inspection of select Cluster Plots aids in identifying problematic SNPs and enables you to manually change calls.

As shown in Figure 114, use the **Type** option above the plot to switch between **Signal** view (left plot), **Log2 Signal** view (not shown), and **Contrast** view (right plot).

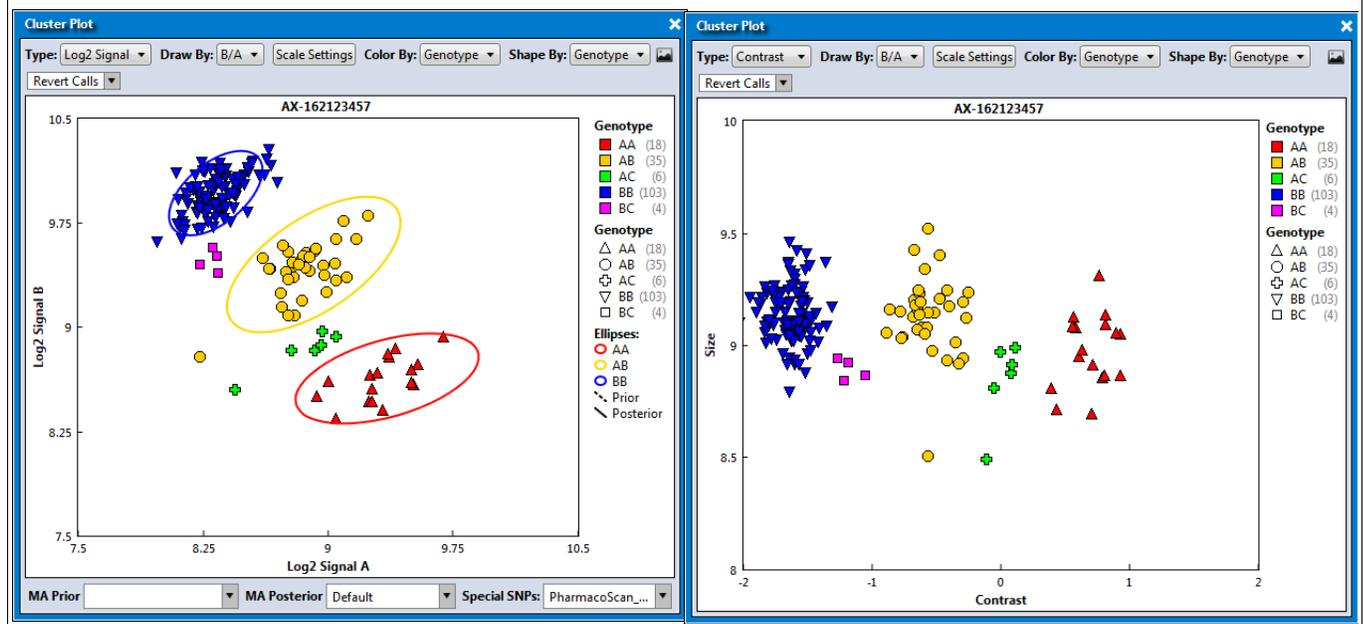
Note: For bi-allele probesets (those that measure two alleles), only the **Contrast** view displays the ellipses that illustrate the prior and posterior knowledge of genotype cluster positions.

Figure 114 Cluster window tab for bi-allelic probesets



Multi-allele probesets measure more than two alleles. Unlike bi-allele probesets, multi-allele probesets are genotyped in the Log2 Signal space. For multi-allele probesets, the ellipses that illustrate the prior and posterior knowledge of genotype cluster positions are only displayed when selecting plot Type **Log2 Signal**. See Figure 115 on page 90 for plot examples of a multi-allele probeset.

Figure 115 Cluster window tab for a multi-allele probeset

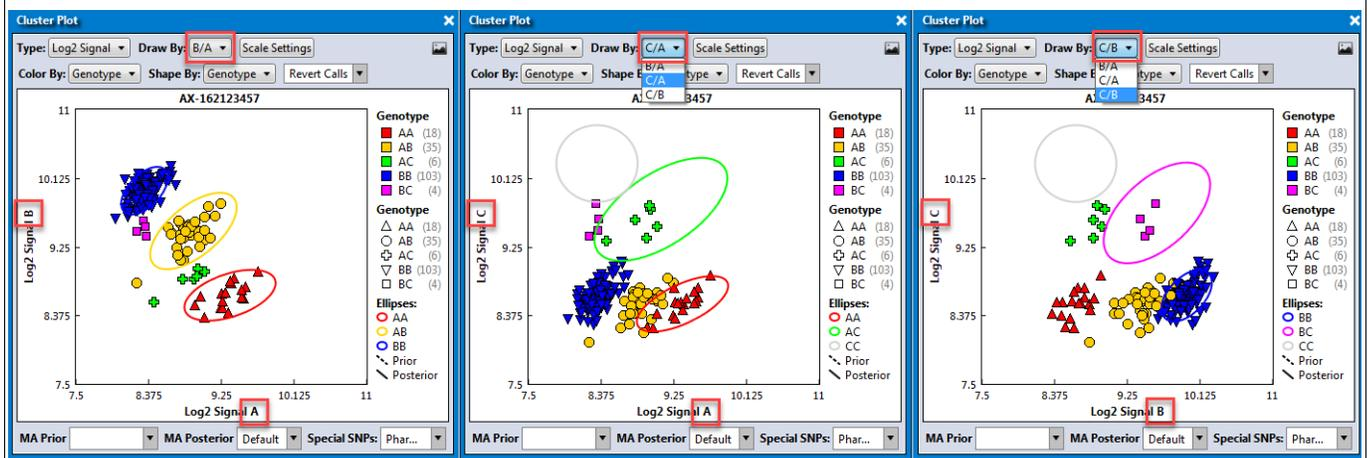


Note: Multi-allele probesets in the SNP Summary Table can be more easily identified by adding the annotation field Allele_Count and/or Ordered_Alleles. If a probeset has Allele_Count of 3 or greater, it is a multi-allele probeset.

Multi-allele probesets measure more than two alleles, and so more than two allele signal channels are present. The cluster plot only shows data for two allele signal channels. If multi-allele probesets are present in the current results, then the top of the cluster plot will also display the Draw By menu.

See Figure 116 for an example of using Draw By to select different pairs of allele signal channels for plotting. The default Draw By option is selected based on the probeset's most common genotype allele calls in the data set.

Figure 116 Cluster window tab - Using Draw By to select different pairs of allele signal channels for plotting.



Using the Cluster Plot

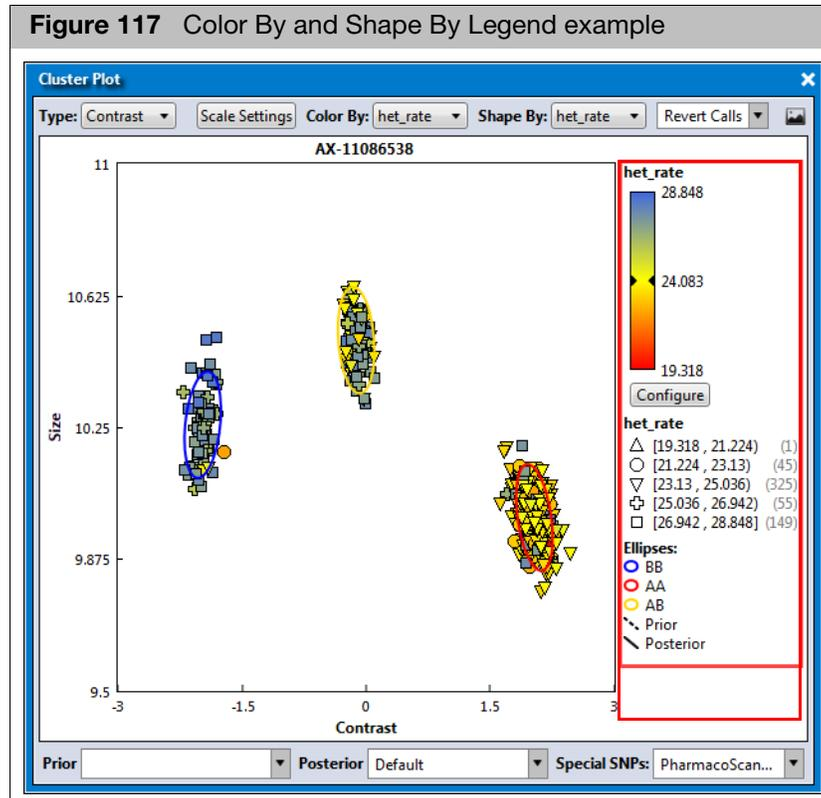
Displaying a SNP Cluster Plot that Corresponds with a SNP

1. In the SNP Summary Table, click on row (SNP) of interest.

Note: Use the arrow keys on the keyboard to toggle through the list. As you toggle through the list, the Cluster Plot auto-updates to match your selected SNP.

2. Click the **Color By** and **Shape By** drop-down menus to select the combination (X and Y axis) view you want. See [Table 12 on page 154](#) for Color By and Shape By definitions. Note: Your imported sample attributes are also available for use within the Color By and Shape By drop-down menus.

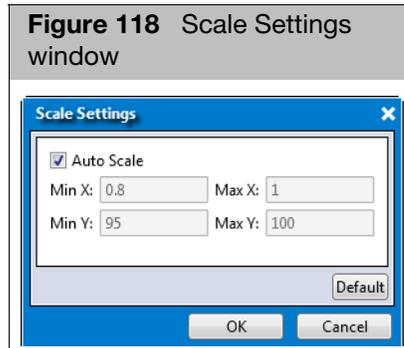
The appropriate legend appears within the plot. (Figure 117)



Setting New Scale Setting Ranges

1. Click **Scale Settings**.

The following window appears. (Figure 118)



By default, the window displays your current range values.

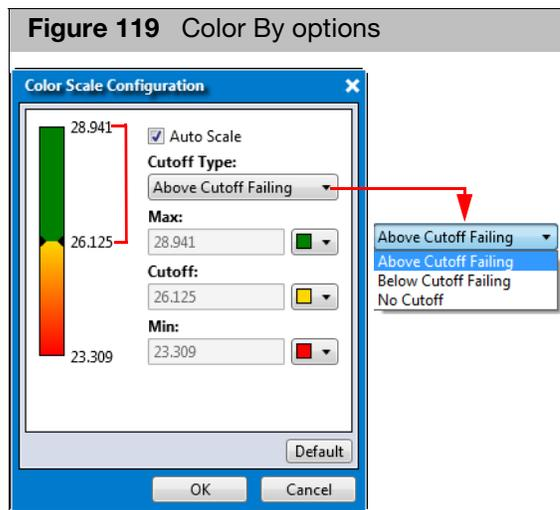
2. Enter your new ranges in the appropriate text fields.
3. Click **OK**.

Click **Default** to return all values back to their factory settings.

Customizing Color By Settings

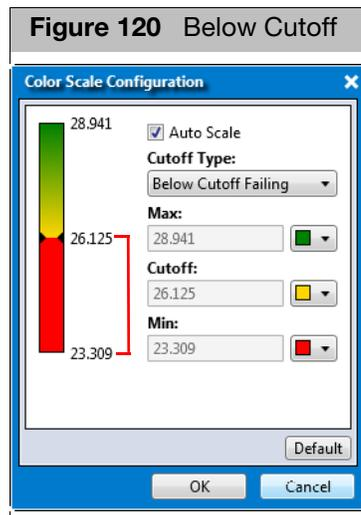
1. Click **Configure**.

The Color Scale Configuration window appears. (Figure 119)

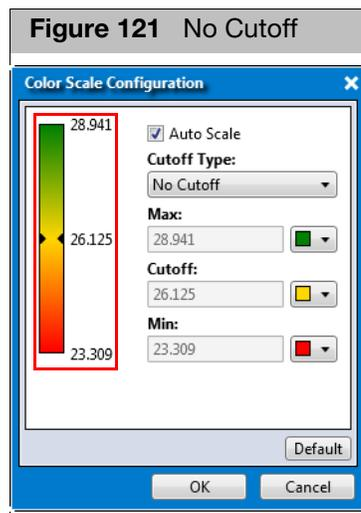


2. Use the provided text fields and color drop-down menus to customize your Color By selection.
 - **Auto Scale** check box (when checked) uses the actual minimum (lower bound) and maximum (upper bound) as your min/max scale. Uncheck the Auto Scale check box to enter your min and max number scales in the provided fields. Note: If the Auto Scale check box remains unchecked, you must enter new scale limits for each subsequent analysis.
 - Click the **Cutoff Type** drop-down menu to select your cutoff preference.
 - **Above Cutoff Failing** - This presents a hard visual cutoff graph of all values that fail ABOVE the Cutoff value entered. The Above Cutoff data is represented by the color defined for Max. (Green in Figure 119)

- **Below Cutoff Failing** - This presents a hard visual cutoff graph of all values that fail BELOW the Cutoff value entered. The Below Cutoff data is represented by the color defined for Min. (**Red** in Figure 120)



- **No Cutoff** - This presents a smooth 3-point gradient of your defined Max, Min, and colors. (Figure 121).



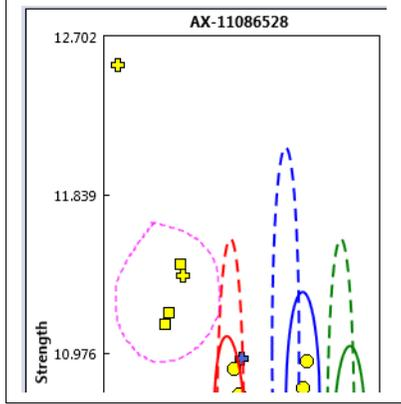
3. Click **OK**.

Your Cutoff preference, entered values, and color selections are now displayed on the graph and saved for future use. If needed, click the **Default** button to revert ALL values back to their factory setting.

Selecting Multiple Samples in a Cluster Plot

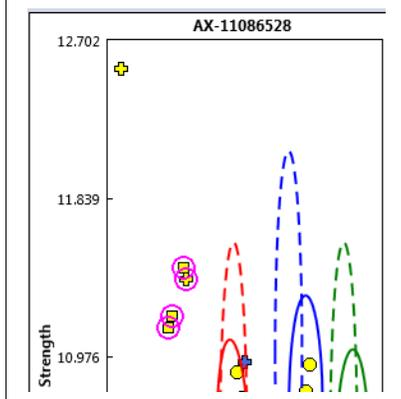
1. Drag the cursor around a group of samples to draw a pink-dotted closed loop around them, as shown in [Figure 122](#).

Figure 122 Selecting multiple samples



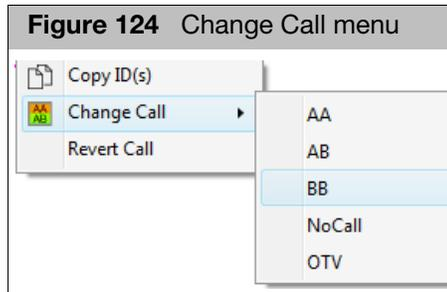
2. Release the mouse button to select the group of lassoed samples. ([Figure 123](#))

Figure 123 Selecting multiple samples



Changing a Sample's Call for a Single SNP

1. Highlight the sample or samples you want to modify, then right-click on them. A menu appears.
2. Click **Change Call**, then move your cursor to the right, then click to select a different call, an OTV (Off Target Variant), or No Call. (Figure 124) Note: If the array supports CN-aware Genotyping, additional Call Code choices appear.



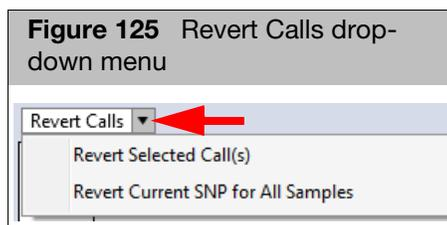
The Call is now changed, but not the position. The image may or may not change, as it depends on the Color By and Shape By options you selected.

Reverting a Single Call

1. Single-click to highlight the Call you want to revert back, then right-click on it. A menu appears.
2. Click **Revert Call**.

Reverting Multiple Calls

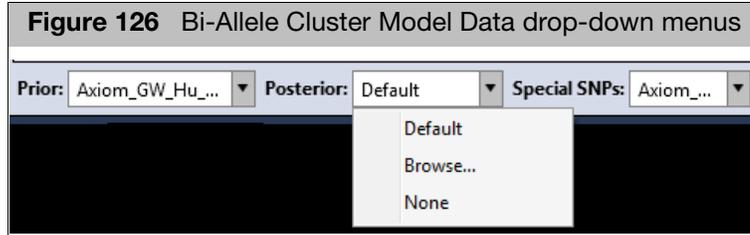
1. Drag the cursor around a group of samples to draw a pink-dotted closed lasso shape around them.
2. Release the mouse button to highlight your selected samples.
3. Click the **Revert Call** drop-down menu (Figure 125), then click to select **Revert Selected Call(s)**.



Displaying Cluster Model Data

By default, the **Prior**, **Posterior**, and **Special SNPs** drop-down selections are preset to best suit the currently displayed Cluster Plot.

1. Click the appropriate drop-down menu (Figure 126), then click to select a new setting.



- If you select **Browse**, an Explorer window appears. Navigate to your folder location as you normally would, then click **Open** to display your data within the graph.
- Selecting **None**, conceals (hides) the selected graph data.

Saving the Current Cluster Plot View

1. Click the **Save Image**  button.
An Explorer window appears.
2. Navigate to where you want to save the .PNG file, enter a filename, then click **OK**.



Allele Translation

About Translations

For supported array types (e.g. PharmacoScan), Axiom Analysis Suite will provide the option to convert (translate) the genotype calls of an important subset of SNPs to functional allele calls using standardized nomenclature wherever possible. The software enables you to:

- Quickly identify possible rare alleles or missing data.
- Identify haplotype and SNP-level sequence variation in the test samples.
- Annotate the reported genotypes across translated SNPs to indicate genomic, mRNA, or peptide changes resulting from any observed variation.
- Predict general gene activity based on detected diplotypes.

See [Appendix B, "About Allele Translation" on page 139](#) for more information.

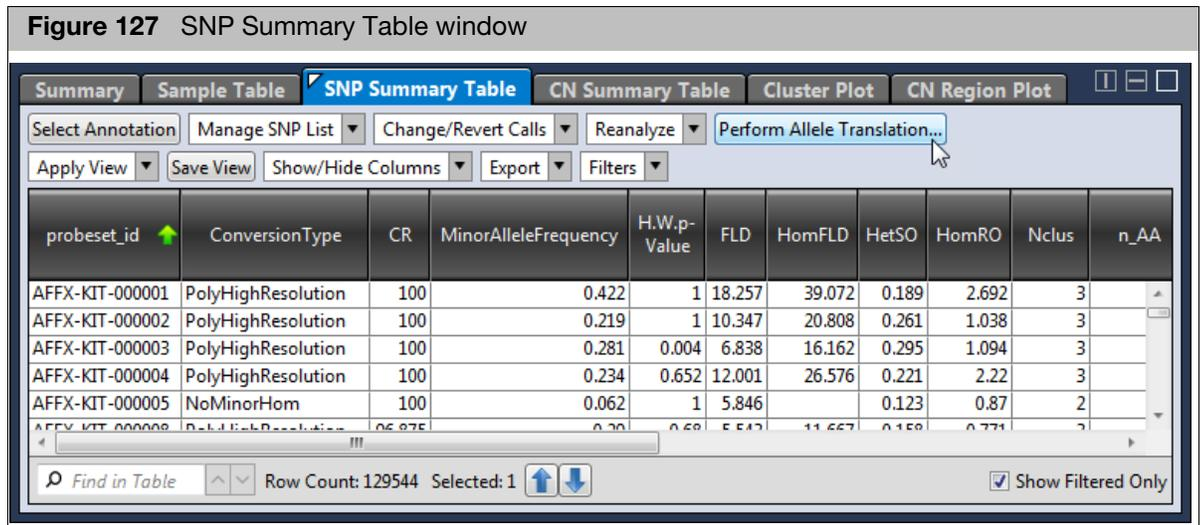
Performing Allele Translation

The Perform Allele Translation operation is available only if the following conditions are met:

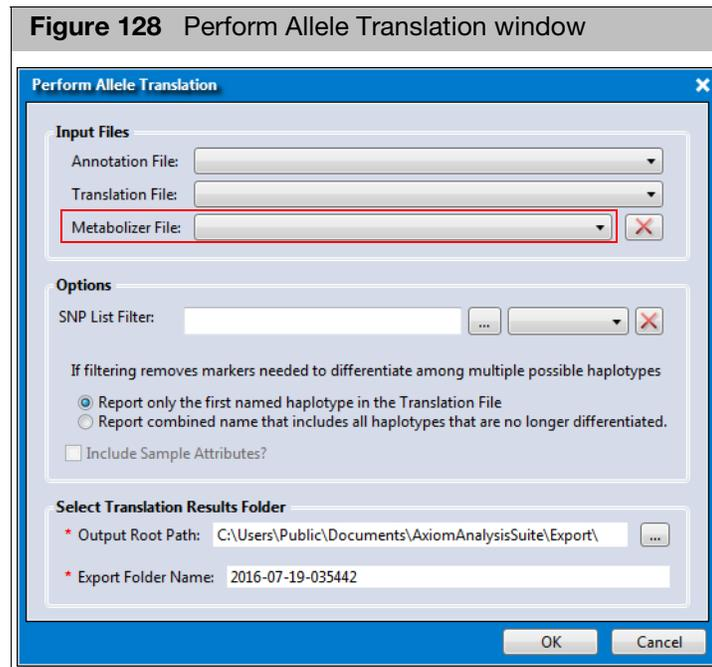
- The library folder must include the same library package (Array Type with revision number) used to generate the batch results.
- Inside the matched library package folder, the files with the extensions ***.dc_annot.csv** and ***.translation** must exist.

To perform an allele translation:

1. From an open batch in the Viewer, navigate to the SNP Summary Table, then click the **Perform Allele Translation** button, as shown in Figure 127.



The Perform Allele Translation window appears. (Figure 128)



2. At the Perform Allele Translation window, use the provided drop-down menus to select an **Annotation File**, **Translation File**, and an optional **Metabolizer File**.
Note: If you disagree with the phenotype interpretations, you may want to leave the Metabolizer File option blank. (Figure 128) A Phenotype report will not be created if this option is not used.

Allele Translation Options

Refer to [Table 1](#) and [Table 2](#) for descriptions of the available **Allele Translation** options.

Table 1 Allele Translation options

Select Options	Description
SNP List Filter	<p>Choose this option to translate only the genotypes of SNPs in a user-specified probeset list. Click the Browse button to select the marker list, or select from SNP lists you're already imported to the batch results.</p> <p>WARNING: If you supplied a custom optional SNP List File for Genotyping in Analysis Settings when setting up your analysis, you should either filter on the same marker list for Allele Translation, or filter on an even more restrictive list. Otherwise, you may see SNPs with NotAvailable calls in your translation reports. The presence of NotAvailable calls may increase the number of reported haplotype possibilities.</p>
Report only the first named haplotype in the translation file (Default)	<p>This option is only relevant IF you filter by a SNP List AND if the SNP List contains some (but not all) of the available probesets in a gene that is allele translated. If both conditions are met, then it is possible that your SNP List excludes a probeset needed to differentiate among two or more named haplotypes in the translation library file.</p> <p>Selecting this option means that only the first haplotype will be reported from the set of possible haplotypes that are non-distinguishable due to probeset exclusion. The haplotypes are ordered by name from left to right in the translation library file.</p> <p>For example, in gene CYP1A1, the *2C haplotype may be differentiated from the *1 haplotype by a variation in a single probeset. If this probeset is omitted, and the data indicates that both *1 and *2C are possible (due to a NoCall at that probeset), then only *1 is reported as a possibility (since *1 is listed before *2C in the translation library file).</p> <p>The disadvantage of selecting this option is that you may be excluding the actual haplotype for a tested sample. The advantage of selecting this option is that you may want to exclude haplotypes that are differentiable only by probesets you have decided not to translate, AND you agree that the selection of which haplotype to report is correct.</p>

Table 1 Allele Translation options

Select Options	Description
Report combined name that includes all haplotypes that are no longer differentiated	<p>This option is only relevant IF you filter by a SNP List AND if the SNP List contains some (but not all) of the available probesets in a gene that is allele translated. If both conditions are met, then it is possible that your SNP List excludes a probeset needed to differentiate among two or more named haplotypes in the translation library file.</p> <p>Selecting this option means that a combined haplotype name will be reported using the set of possible haplotypes that are non-distinguishable due to probeset exclusion.</p> <p>For example, in gene CYP1A1, the *2C haplotype may be differentiated from the *1 haplotype by a variation in a single probeset. If this probeset is omitted, and the data indicates that both *1 and *2C are possible (due to a NoCall at that probeset), then "**1_or_*2C" is reported as a possibility.</p> <p>The advantage of selecting this option is that you are not excluding possible haplotypes. The disadvantage of selecting this option is that the report will include haplotypes that require a variant allele of a probeset you have decided to exclude for translation.</p> <p>Note: This option is only available if you do not need a phenotype report, as phenotyping requires haplotype names to not change (depending on the set of probesets used for translation). To enable this option that excludes the generation of a phenotype report, you must deselect the usage of the metabolizer library file. To do this, click  (right of the Metabolizer File option).</p>
Include Sample Attributes	Click this check box to include sample attributes in the translation reports. This option is enabled if you have imported sample attributes to your analysis results.

Table 2 More Allele Translation options

Select Translation Results Folder	Description
Output Root Path	The path to the output folder for the translation reports and log. Click its Browse button to set the path.
Export Folder Name	The new folder name for the translation results, whose default name has the date-time format YYYY-MM-DD_HHMMSS_translations. Confirm the default or enter a new folder name.

- After completing the Allele Translation selections, click **OK**.
When the translation is finished, an Explorer window appears displaying the folder containing the translation reports and corresponding run log.

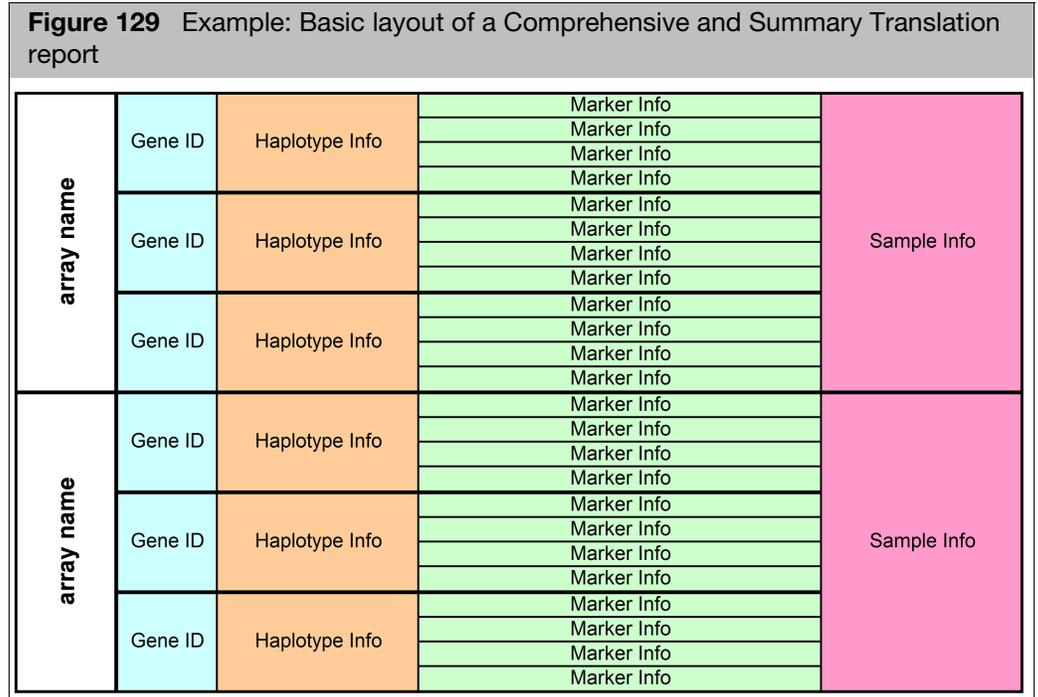
Translation Reports

The allele translation operation creates the following reports and supporting files:

- **Comprehensive Translation** - Displays one row per translated SNP for each sample. Provides information on each SNP in addition to haplotype calls.
- **Summary Translation** - An abbreviated version of the Comprehensive report, which displays at least one row for every translated gene for each sample. It also includes rows for every genotype where the translation identifies a variant call. It also includes rows listing SNPs with missing data. In the Summary report only, if no SNPs responsible for functional changes report a variant allele, then information for those SNPs is replaced with a comment to this effect. If a copy number state of zero is indicated, then information for SNPs in that gene is replaced with a comment to this effect, and the copy number haplotype code is reported in the Known Call field.
- **Phenotype Translation** - Displays one row per phenotyped gene for each sample, based on the diplotypes from the source Comprehensive Translation report
- **Uncalled probeset list** - A list of probesets with NoCall genotype calls from SNPs used for translation. This probeset list can then be imported into the Batch, so that you can filter the SNP Summary Table by the uncalled.ps SNP list. This provides a quick way to review the cluster plots of probesets with missing data, and possibly edit the calls directly to "fill in" the missing data in preparation for another round of allele translation reports.
- **Uncalled** - A list of NoCall genotype calls from SNPs used for translation. This report is useful for identifying samples and SNPs for follow up genotyping.
- **MD5** - An electronic signature that can be used to verify that the comprehensive and phenotype reports have not been modified. Interested users can contact devnet@affymetrix.com for information on accessing tools to verify the integrity of the translation results files.
- **Log file** - A list of messages generated by the software as the data is processed. This file is useful for troubleshooting errors.

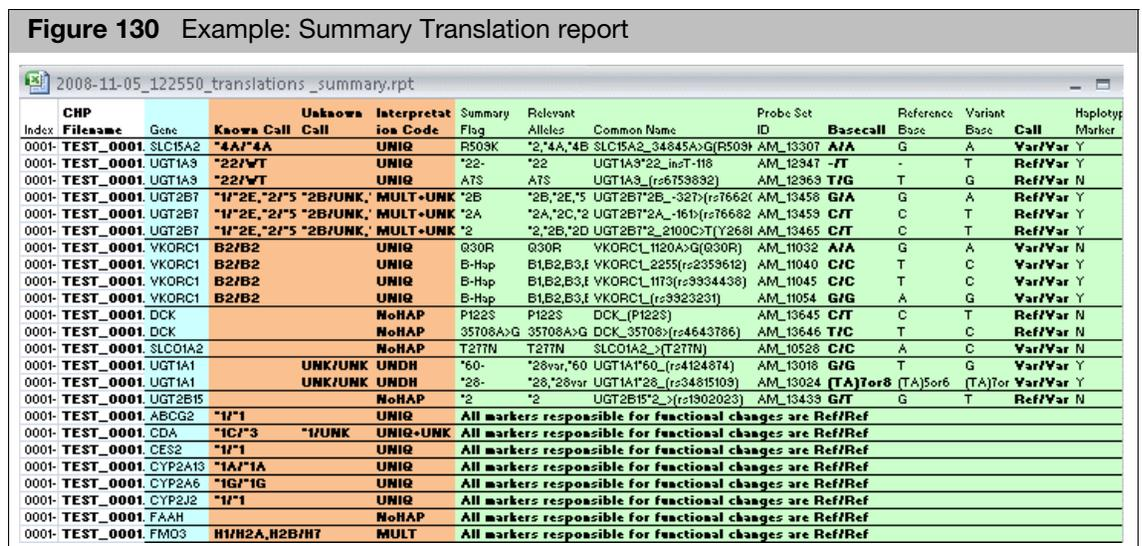
Comprehensive and Summary Translation Report

The basic layout of this report is shown in Figure 129.



Summary Translation Report

To make this report easier to read, bold and regular fonts are used. For example, genotype calls are represented in bold, as shown in Figure 130.



Phenotype Translation Report

The basic layout of this report is shown in Figure 131.

Figure 131 Example: Phenotype Translation report

array name	Gene ID	Phenotype Info	Haplotype Info	Sample Info
	Gene ID	Phenotype Info	Haplotype Info	
	Gene ID	Phenotype Info	Haplotype Info	
	Gene ID	Phenotype Info	Haplotype Info	
array name	Gene ID	Phenotype Info	Haplotype Info	Sample Info
	Gene ID	Phenotype Info	Haplotype Info	
	Gene ID	Phenotype Info	Haplotype Info	
	Gene ID	Phenotype Info	Haplotype Info	

Phenotype Report

The basic layout of this report is shown in Figure 132.

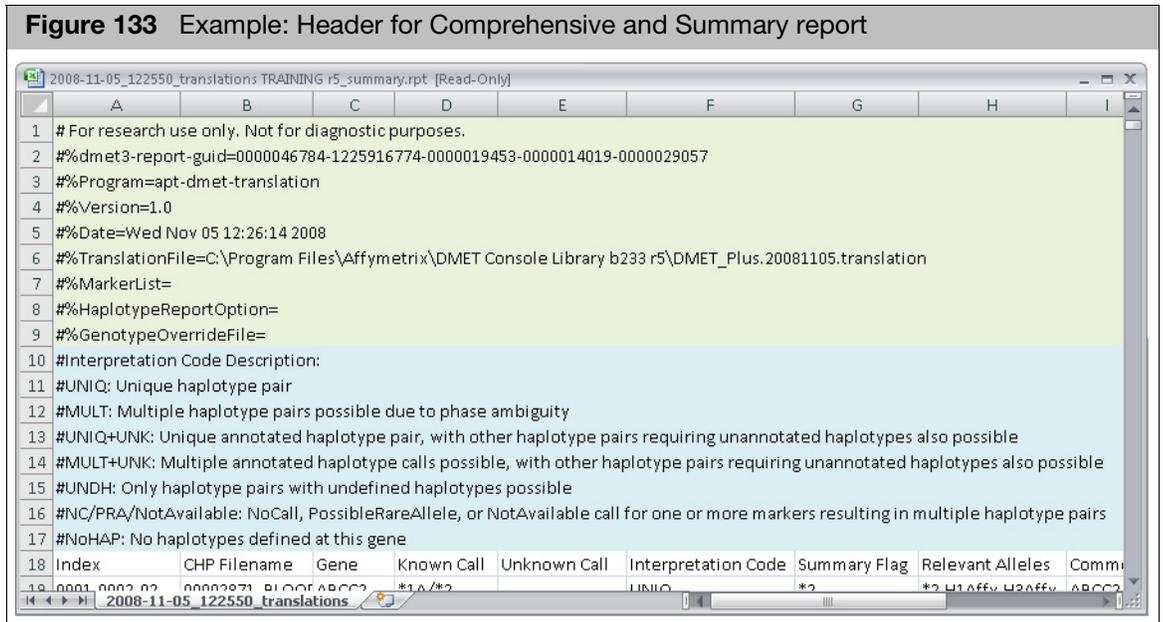
Figure 132 Example: Phenotype report

Index	CHP File	Gene	Phenotype Call	Gene Activity	Known Call	Unknown Call	Interpretation Code
0001-0020	test_01.c	CYP1A2	EM	normal/normal	*1F/*1F		UNIQ
0001-0022	test_01.c	CYP2A6	EM	normal/normal	*1/*1		UNIQ
0001-0024	test_01.c	CYP2B6	EM	normal/normal	*1/*1		UNIQ
0001-0029	test_01.c	CYP2D6	PM	none/none	*5/*5		UNIQ
0002-0020	test_02.c	CYP1A2	EM_or_IM	normal/reduced	*1A/*1L,*1C/*1F		MULT
0002-0022	test_02.c	CYP2A6	EM_or_IM	normal/reduced	*1/*17		UNIQ
0002-0024	test_02.c	CYP2B6	EM_or_IM	normal/reduced	*1/*6	*4/UNK	UNIQ+UNK
0002-0029	test_02.c	CYP2D6	EM_or_IM	normal/reduced	*2/*29	*2/UNK,*29/UNK	NC/PRA/NA
0004-0020	test_04.c	CYP1A2	EM	normal/normal	*1A/*1A		UNIQ
0004-0022	test_04.c	CYP2A6	EM	normal/normal	*1/*1		UNIQ
0004-0024	test_04.c	CYP2B6	IM	normal/none	*1/*18		UNIQ
0004-0029	test_04.c	CYP2D6	EM	normal/normal	*1/*1		UNIQ
0005-0020	test_05.c	CYP1A2	EM	normal/normal	*1A/*1F		UNIQ
0005-0022	test_05.c	CYP2A6	EM	normal/normal	*1/*1		UNIQ
0005-0024	test_05.c	CYP2B6	IM	reduced/reduced	*6/*6		UNIQ
0005-0029	test_05.c	CYP2D6	IM	normal/none	*2/*4	*1/UNK,*2/UNK,*4/UNK	NC/PRA/NA
0009-0020	test_09.c	CYP1A2	EM	normal/normal	*1F/*1F		UNIQ
0009-0022	test_09.c	CYP2A6	EM	normal/normal	*1/*1		UNIQ
0009-0024	test_09.c	CYP2B6	IM	reduced/reduced	*6/*6		UNIQ
0009-0029	test_09.c	CYP2D6	IM	normal/none	*2/*4	*1/UNK,*10/UNK,UN	UNIQ+UNK

Opening Translation Report in MS Excel

1. Use Windows Explorer as you normally would to navigate to the export folder with the translation results.
2. Double-click the report (.rpt) to be viewed. You may be asked choose an application to open the report. Select Microsoft Excel, then step through the Text Import Wizard (use the tab-delimited default options).

The report header includes basic information that helps track study data and definitions of interpretation codes, as shown in [Figure 133](#).



Available Report Fields and Descriptions

Array Tracking

Refer to [Table 3](#) for descriptions of the available **Array Tracking** fields.

Table 3 Array Tracking fields

Array Tracking fields	Description
Index	A row index in the format: [filename index]-[gene index within filename]-[Probe Set ID index within gene]. This field can be parsed for sorting or row filtering. For the phenotype report, the index is shortened to [filename index]-[gene index within filename]
Filename	Name of the sample file.

Gene-specific

Refer to [Table 4](#) for descriptions of the available **Gene-specific** fields.

✗ IMPORTANT! Haplotypes are not reported for genes whose Interpretation Code is **NoHap**. The fields described in the table below will therefore be empty for these genes. The exception is if the gene reports a gene deletion, in which case the associated haplotype names are reported.

Table 4 Gene-specific fields

Gene-specific fields	Description
Associated Gene	Gene symbol
Phenotype Call	<p>In the Phenotype report, the predicted phenotype given the supplied Known Call diplotypes. Multiple comma-separated phenotypes are reported when multiple Known Call diplotypes are associated with different phenotypes. Most genes use the following terminology when the default metabolizer library file is selected:</p> <ul style="list-style-type: none"> • UM = ultra-rapid metabolizer • RM = rapid metabolizer • NM = normal metabolizer • IM = intermediate metabolizer • PM= poor metabolizer <p>Variations on these terms also exist to describe some level of uncertainty:</p> <ul style="list-style-type: none"> • NM_or_IM = normal or intermediate metabolizer • IM_or_PM =intermediate or poor metabolizer • Not_PM = not a poor metabolizer • unknown = unknown metabolizer state <p>Some genes use different phenotype terms to be consistent with literature usage. Refer to the header of the phenotype report for additional information.</p> <p>Users are responsible for reviewing the *.metabolizer library file for accuracy! Users may modify the *.metabolizer file as needed, and are not restricted to this terminology. Refer to "Diplotype to Phenotype Translation" on page 144 for more information.</p>

Table 4 Gene-specific fields

Gene-specific fields	Description
Gene Activity	<p>In the Phenotype report, the predicted pair of gene activities given the supplied Known Call diplotypes. Multiple comma-separated activity pairs are reported when multiple Known Call diplotypes are associated with different activity pairs. Most genes use the following terminology when the default metabolizer library file is selected:</p> <ul style="list-style-type: none"> • increased = increased gene function • normal = normal gene function • reduced: reduced gene function • no = no gene function • unknown = unknown or uncertain gene function <p>Some genes use different phenotype terms to be consistent with literature usage.</p> <p>Users are responsible for reviewing the *.metabolizer library file for accuracy! Users may modify the *.metabolizer file as needed, and are not restricted to this terminology. Refer to "Diplotype to Phenotype Translation" on page 144 for more information.</p>
Known Call	<p>Haplotype pairs (diplotypes) identified in the gene of interest. When more than one pair of haplotypes is implicated (due to phase ambiguity in compound heterozygous samples), the reported diplotypes are separated by a comma.</p>

Table 4 Gene-specific fields

Gene-specific fields	Description
Unknown Call	<p>When the gene table includes haplotyping SNPs and a complete diplotype pair cannot be identified in a sample, one or more unknown haplotypes is assumed. This is designated as UNK in the report. Multiple haplotype pairs (diplotypes) that have unknown alleles are separated with a comma in this field. An example record might be in the format: *2/UNK,*13/UNK,*24/UNK,*32/UNK to indicate that there are at least 4 defined alleles consistent with the data, but each would require matching to a haplotype pattern that does not exist in the translation library file.</p>
Interpretation Code	<p>This diplotype interpretation code indicates whether one and only one unique haplotype pair is consistent with the data (UNIQ), whether there are multiple haplotype pairs consistent with the observed genotypes (MULT) and whether these are observed in conjunction with other unknown haplotypes (UNIQ+UNK or MULT+UNK). Additional codes indicate that no known haplotype pairs have been identified (UNDH) or if there is missing data leading to additional haplotype possibilities. The missing data could be NoCall or NotAvailable (NC/PRA/NA). The PossibleRareAllele call is not used by Axiom Analysis Suite.</p> <p>For genes for which copy number state is available, the following interpretation codes may also appear:</p> <ul style="list-style-type: none"> • CN_HybridLoss = Partial gene deletion is detected, so haplotype pair calling is not available. • CN_HybridGain = Partial gene duplication is detected. If a non-wild-type allele is detected, the software can't determine whether the variant allele is on the partial copy of the gene. • CN_Gain = Gene duplication is detected, but the software cannot determine which haplotype(s) are duplicated. For this reason you won't see xN nomenclature in the reported haplotype pairs. • CN_NoCall = Copy Number state not reported, so there is less confidence in the reported haplotype pairs. • CN_Error = Genotypes from multiple Copy Number states are detected, so haplotype pair calling is not available. This can happen if you manually edit a genotype call in such a way as to change the SNP's copy number state. For example, if you change a "NoCall_1" genotype (a NoCall for a CN=1 sample) to a "BB" CN=2 genotype, its copy number has changed. This would cause allele translation problems if other SNPs for this sample for the same gene have genotypes of a different copy number state.

Marker-specific

Refer to [Table 5](#) for descriptions of the available **Marker-specific** fields.

✗ IMPORTANT! Marker annotations will always be the same throughout the reports for a given SNP. Fields in the Marker Information section of the Comprehensive and Summary reports include the biological information at the SNP level, along with the interpreted genotypes identified in each sample

Table 5 Marker-specific fields

Marker-specific fields	Description
Summary Flag (marker annotation)	<p>This annotation field contains an abbreviated name when structural or functional differences are known to result with mutations at the SNP locus. For triallelic SNPs, there may be more than one flag. For example, ABCB1_c.2677G>T>A(A893S or T) marker is triallelic and can result in different function changes in the protein. The two summary flags for the marker are thus reported "A893S,A893T".</p> <p>This flag is N (No) if the marker is not defining for a haplotype, and its genotype doesn't affect the structure or expression of the gene product. If the flag is N, the marker will not appear in the Summary report unless it reports NoCall.</p>
Relevant Alleles (marker annotation)	<p>This annotation field is the full listing of haplotype-based alleles defined in the gene table that contain the variant version of the marker. For non-haplotype-based SNPs, this is an abbreviated name indicating the protein change that results when the variant base is present.</p>
Common Name (marker annotation)	<p>A SNP identifier describing either the gene location, coding change or dbSNP rsID for the SNP. The Common Name is retrieved from the translation library file, and may not be the same as the Common Name seen within Axiom Analysis Suite (which instead uses the Common Name in the *.annot.db library file)</p>
Probe Set ID	<p>Unique identifier for the SNP.</p>
Basecall	<p>The observed bases, also known as the "raw" genotypes.</p>
Reference Base	<p>This field generally indicates the more common allele in biallelic SNPs. Certain genes use a particular GenBank entry as the "Reference genome" and the observed allele at each marker across the gene is then reported as Reference.</p>
Variant Base	<p>These are the alternate alleles for each SNP. When there is more than one variant allele (e.g. triallelic SNPs) the alternate alleles are reported together and separated by a comma (e.g. A,T).</p>
Call	<p>The first level of translation of the Basecall field, replacing the individual nucleotide calls with the associated reference (Ref) or variant (Var) allele state. For Basecalls associated with copy numbers less than 2, this field will show haplotype names as needed. This field will be empty if the Basecall value has an unclear call.</p>
Haplotype Marker (marker annotation)	<p>Differentiates SNPs used to make haplotype calls or single-marker variant calls.</p> <ul style="list-style-type: none"> • Y = A flag to indicate that the Allele translation algorithm will match allele variants in blocks of SNPs defining haplotypes in the gene tables. Called haplotypes are reported in the "Known Calls" and Unknown Calls of the report. • N = A flag to indicate that the haplotype background of a variant is not known. Genotyping results for these SNPs are only reported in the "Call" column.
Change for Variant (marker annotation)	<p>Amino acid substitution or other structural change (such as splicing variant, promoter mutation, Frame shift mutation, etc.) caused by the presence of the variant allele.</p>

Table 5 Marker-specific fields

Marker-specific fields	Description
cDNA Change (marker annotation)	Location of the mutation on a reference mRNA sequence.
Genome Position (marker annotation)	This is the chromosomal position of the mutation.
dbSNP RS ID (marker annotation)	The dbSNP identifier for the marker.

Tracking Edited Genotype Calls

Refer to [Table 6](#) for descriptions of the available fields for tracking edited genotype calls.

X IMPORTANT! Fields for tracking genotyping changes are recorded in the Change Tracking portion of the translation reports (located immediately before the User Defined Sample Information section of the report).

Table 6 Fields for tracking edited genotype calls

Tracking Edited Genotype Calls	Description
Original Basecall	This field reports the original genotype for the sample. The field is not empty when the user has edited the call within Axiom Analysis Suite.
Override Comment	This field reports edited in AxAS if the call has been edited within Axiom Analysis Suite.

Uncalled Report

Refer to [Table 7](#) for descriptions of the available fields for uncalled reports.

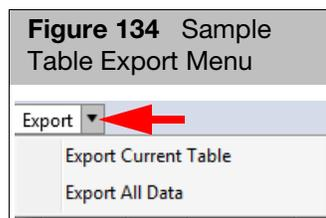
X IMPORTANT! The Reference Allele and Variant Allele values of the Uncalled report are the same as the Reference Base and Variant Base values of the Comprehensive and Summary reports.

Table 7 Uncalled Report

Uncalled Report	Description
Filename	Name of the sample file.
Gene	Gene symbol.
Common Name	The Common Name defines the gene and positional information about the genetic change tested with the probeset.
Basecall	The Uncalled report contains all the NoCall genotypes from SNPs used for allele translation.
Override Comment	User-specified annotation field that enables an audit trail of the source of genotyping results done outside of Axiom Analysis Suite. In the Uncalled report, this field is primarily useful when creating a copy of this file to use as an Override file. The Override file may be useful for Affymetrix Power Tool users.
Reference Allele	Reference base indicates the allele in a reference sequence known to be present at this genetic location. Generally this is the more common allele at SNPs with low minor allele frequency (<1%).
Variant Allele	The variant base(s) defined by the marker are alternative known genotypes known to be present at this genetic location. For triallelic SNPs, the reporting format is: A,T for ABCB1_68883G>T(S893A) because two specific mutations are known to occur at this genomic location (G>A and G>T).

Using the Sample Table Export Options

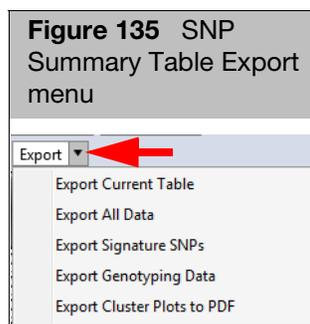
1. Click the **Export** drop-down.
Your Export options appear. (Figure 134)



2. Click **Export Current Table** or **Export All Data**.
A **Save As** window appears.
3. Click on an existing folder or click **New Folder** to choose a new save location.
4. Type a filename for the table, then click **Save**.
The table data is now saved as a tab-delimited .txt file.

Using the SNP Summary Table Export Options

1. Click the **Export** drop-down.
The following window appears: (Figure 135)



Exporting the Current Table

To export the columns and rows currently displayed in the table:

1. Click **Export Current Table**.

An Explorer window appears.

2. Enter a filename, then click **Save**.

The current table data is now saved and exported as a tab-delimited text file.

Exporting All Data

To export all columns and rows, including hidden and filtered data:

1. Click **Export All Data**.

An Explorer window appears.

2. Enter a filename, then click **Save**.

All data in the table (displayed or not) is saved and exported as a tab-delimited text file.

Exporting Signature SNPs

Use this option to export only the signature SNPs in your data.

1. Click **Export Signature SNPs**.

An Explorer window appears.

2. Enter a filename, then click **Save**.

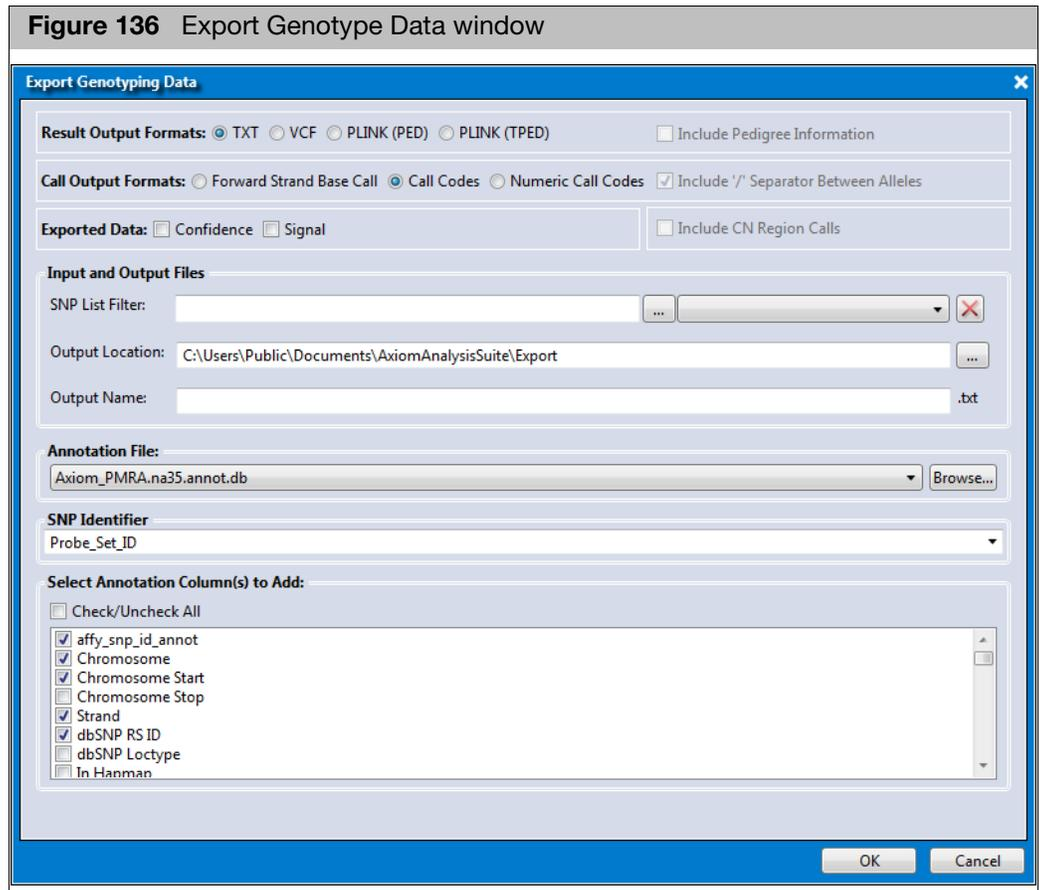
All data is now saved a tab-delimited text file.

Exporting Genotyping Data

 **IMPORTANT!** Not all options are available and are dependent on the export format you select and its applicable format restrictions.

1. Click **Export Genotyping Data**.

The following window appears: [\(Figure 136\)](#)



Result Output Formats

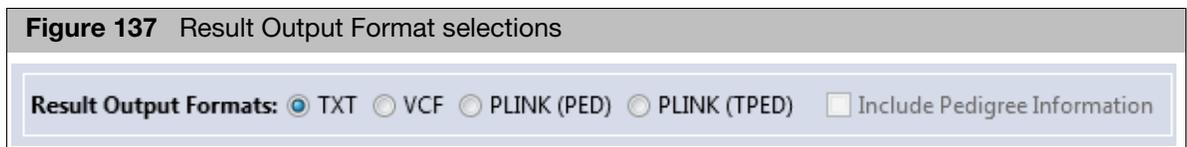


Figure 137 Result Output Format selections

1. Click the radio button to select the Result Output Format you want use.
2. Optional: If you selected a PLINK format, make sure you click the **Include Pedigree Information** check box. Not checking this box may require special handling (within PLINK) to make your exported output work properly.

Note: PLINK format specifies that all markers be biallelic, therefore multi-allele probesets will not be exported to PLINK files.

✗ IMPORTANT! If you click the Include Pedigree Information check box with your PLINK format, make sure your Sample Attributes include an Index ID and Pedigree Information (Family ID, Individual ID, Father ID, Mother ID, Sex, and Affection Status).

3. If you have used OTV Caller, some of the calls may be OTV. To export these OTV calls, click the **TXT** and **Call Codes** radio buttons or **TXT** and **Numeric Call Codes** radio buttons. Note: For all other formats, OTV calls are treated as No Calls.

Call Output Formats

1. Click the radio button to select the appropriate Call Output Format (Figure 138) you want to use.



If you select Forward Strand Base Call, you have the option of not including the '/' separator between alleles. Note: Excluding separators is NOT advised when exporting probesets with multi-base alleles.

✗ IMPORTANT! If exporting Numeric Call Codes, the exported data file's header rows includes a table mapping numeric call code to call code. The format of these header rows is:

```
#%%call-code-N=call_code:numeric_call_code:code_ploidy
```

The assignment of numeric call codes to call codes is NOT guaranteed to be the same across all supported array types.

Exported Data Selections

1. Click inside the check box(es) to check the additional type(s) of Exported Data (Figure 139) you want to include.

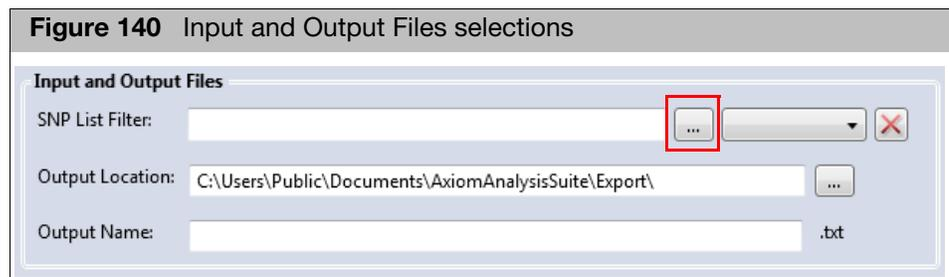


Note: Multi-allele probesets will not export signal values. Normalized signal values for multi-allele and bi-allele probesets are available in the AxiomGT1.summary.txt file inside the batch results folder.

If you selected the VCF output format and copy number analysis was performed, you have the option to not include CN Region Calls from the VCF file.

Input and Output Files

(Figure 140)



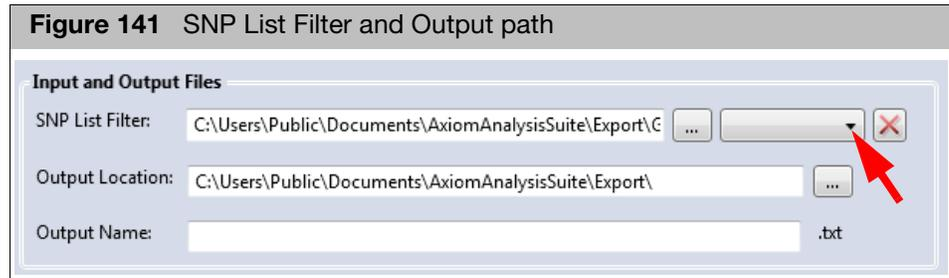
SNP List Filter (Optional)

1. To restrict the output to a list of SNPs (probeset_IDs) contained in a file, click the SNP List Filter field's **Browse** button (Figure 140).

An Explorer window appears.

2. Navigate to the SNP Filter List location, then click **Open**.

The SNP Filter List field is now populated. (Figure 141)



3. Click the SNP List Filter's drop-down menu to view and select a previously saved SNP List. (Figure 141)

Output Location (Required)

1. Click the Output Location field's **Browse** button.

An Explorer window appears.

2. Navigate to an output location, create a new folder if needed, then click **Select Folder** button.

The Output Location path is displayed. (Figure 141)

Output Name (Required)

1. Use the output name already in the Output Name field, or click inside the field to enter a new name. Note: Your output name's file extension reflects the Results Output Format you selected in Step 1.

Changing the SNP Identifier

✗ IMPORTANT! Only SNPs that have a value for the selected annotation are exported. For example, markers that have a dbSNP RSID are exported, while markers without a dbSNP RSID are not exported.

If the selected SNP Identifier has more than 1 probeset mapped to it, it will have multiple entries in the exported file.

1. Click the drop-down arrow, then click to select the SNP Identifier you want to use. (Figure 142)



Changing the Current Annotation File (Optional)

1. To change the currently displayed Annotation File, click the Annotation File field's **Browse** button (Figure 143).

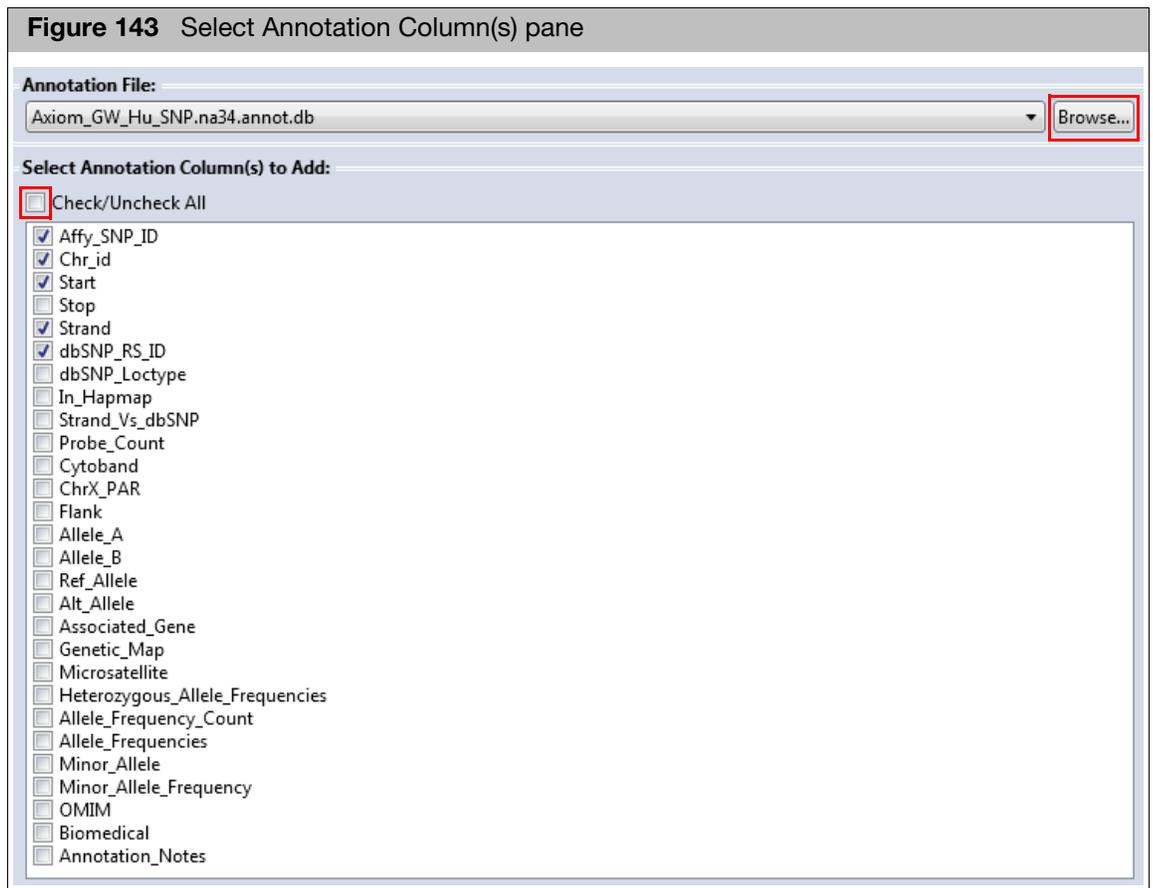
An Explorer window appears.

2. Navigate to the appropriate Annotation File location, then click **Open**.

Your newly selected Annotation file is displayed.

Adding and Removing Annotation Columns

1. Click the check box next to the Annotation Column(s) you want to add to your format results or click to uncheck/remove a column. If you want to add or remove ALL available annotations, click the **Check/Uncheck All** check box. (Figure 143)



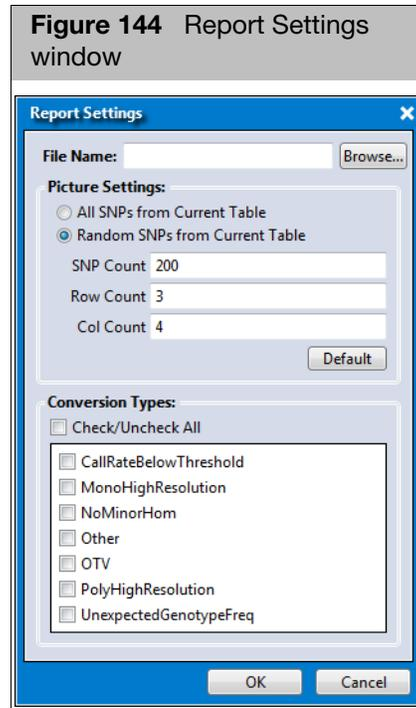
2. After the Export Genotype Data form is complete, click **OK**.
3. Your newly exported data now reside in the output location you defined in [Step 1 on page 115](#).

Exporting Cluster Plots to PDF

This exporting option generates a specific number of cluster plots from the selected Conversion Type choices.

1. Click **Export Cluster Plots to PDF**.

The following window appears: (Figure 144)



2. Click **Browse**.

An Explorer window appears.

3. Navigate to a desired location, then enter a name for your PDF report.
4. Click **Save**.

You are returned to the **Report Settings** window.

5. In the Picture Settings section, click either:

- All SNPs from Current Table
- Random SNPs from Current Table

6. Click inside the applicable **Count** field(s) to enter how many cluster pots you want to export.

Click the **Default** button to revert your modified counts back to their original states.

7. Click to check a Conversion Type(s) you want to add to the report or click the **Check/Uncheck All** check box, then click **OK**.

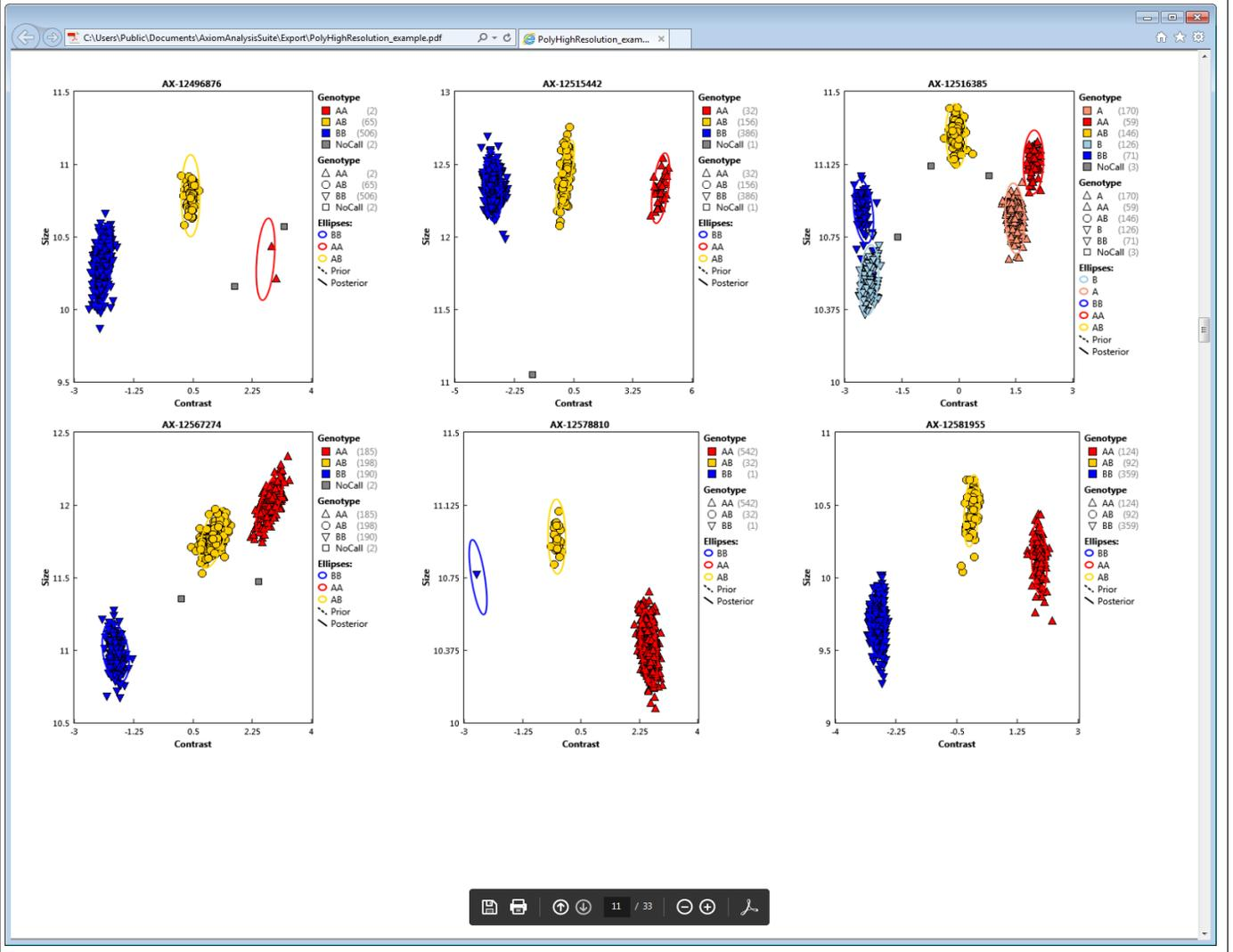
A Please Wait message and progress bar appear. Allow several minutes if multiple Conversion Type PDF report(s) were selected.

An Explorer window (where you saved the PDF Report location in Step 3) appears.

8. Double-click on the PDF Report you want to view.

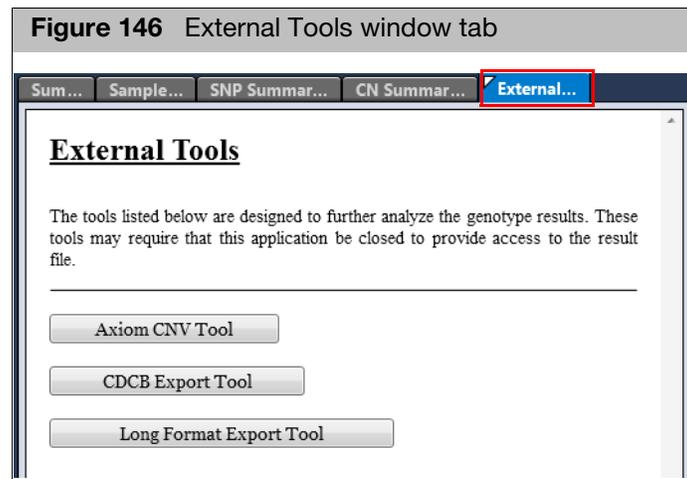
The PDF Report opens. (Figure 145)

Figure 145 PDF Report example



Note: One plot is made for each probeset. For multi-allele probesets, the most informative pair of allele signal channels in the data set is used for plotting, based on the distribution of genotype calls.

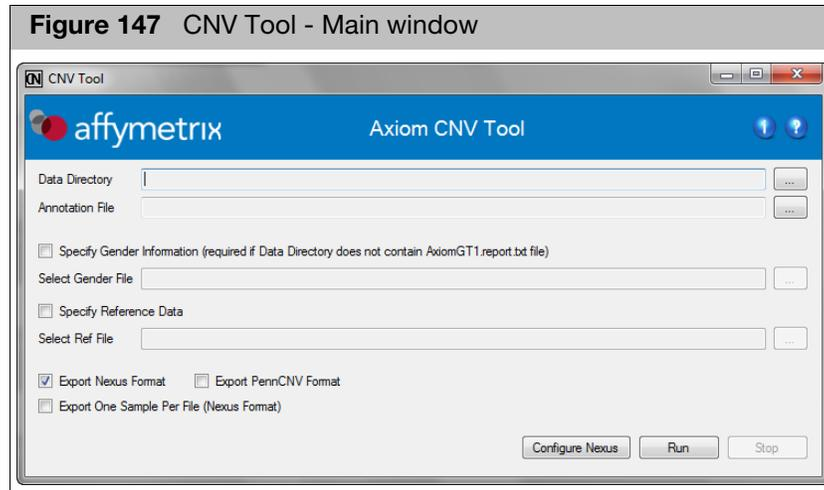
Axiom Analysis Suite auto-detects any previously installed application that may further analyze your genotype results. When a compatible application is detected, an **External Tools** window tab is generated, as shown in [Figure 146](#).



X IMPORTANT! The Axiom Analysis Suite can be left open/active if you are using an external application to further view your data. However, you must exit and close the Axiom Analysis Suite before editing its data with an external application.

Axiom CNV Tool 1.1

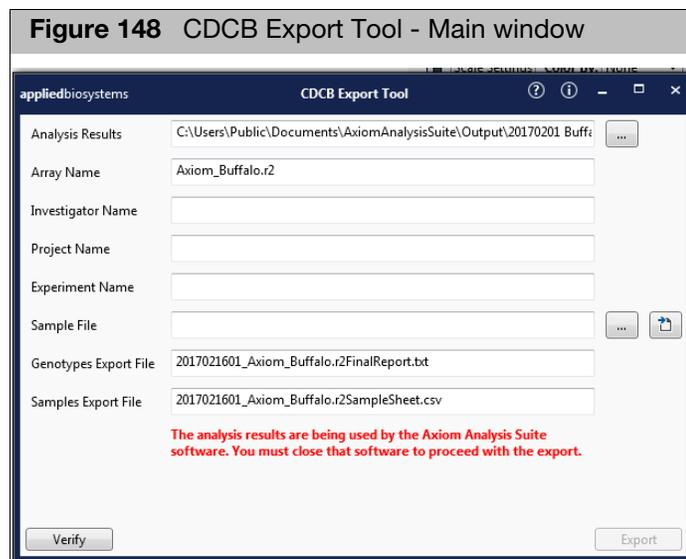
1. Click the **Axiom CNV Tool** button.
The application opens. (Figure 147)



For details on how to use the Axiom CNV Tool 1.1 application, refer to its User Guide (P/N 703216).

Council on Dairy Cattle Breeding (CDCB) Export Tool

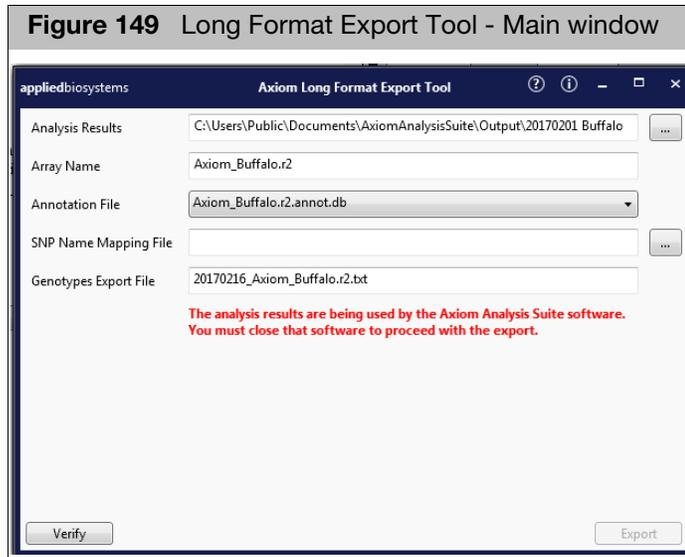
1. Click the **CDCB Export Tool** button.
The application opens. (Figure 148)



For details on how to use the CDCB Export Tool application, refer to its User Guide (P/N 703465).

Axiom Long Format Export Tool

1. Click the **Long Format Export Tool** button.
The application opens. (Figure 149)



For details on how to use the Long Format Export Tool application, refer to its User Guide (P/N 703455).



Predefined Region CN Analysis Variations

Predefined Copy Number Analyses

If the array content and library package supports it, the Genotyping and Best Practices workflows perform predefined region copy number analysis. Copy number analysis is restricted in that the copy number regions are pre-specified, with a fixed set of probesets used to estimate a single copy number state for each region. This form of copy number analysis is done within the Genotyping and Best Practices workflows.

Axiom Analysis Suite supports two variations of restricted copy number analysis. These analysis variations are determined by the array library package and are compared in [Figure 150](#).

Figure 150 Table showing workflow analysis variations that support copy number

Workflow Stage		Analysis Description	Analysis Variation		
			Typical Axiom	Predefined CN Genotyping †	CN-Aware Genotyping ‡
Best Practices	Sample QC	Sample QC	yes	yes	yes
		Identify Copy Number (CN) control samples that pass Sample QC	not done	female samples used as CN plate controls	CN controls identified by comparing measured with reference genotype calls
	Genotyping	Normalize plate signals using CN control samples that pass CN QC. Compute CN state in pre-defined regions. Samples that fail CN QC will report a CN state of NoCall.	not done	yes	yes
		Supply CN results to genotyping engine	not done	no	yes
		Final genotyping	yes	yes, including OffTargetVariant calls	yes, using additional call codes for haploid and zero CN calls
† library package specifies analysis_category = "CN_GT_2" ‡ library package specifies analysis_category = "CN_GT"					

With Predefined CN Genotyping

- CN states will be reported for the predefined regions specified by the library package
- Unless overridden by the user, the control samples used for per-plate signal correction for CN analysis are selected from the set of samples determined to be female and passing sample QC



- Unless overridden by custom library file settings, genotyping probesets will report diploid genotype codes (such as "AA", "AB", "BB", "NoCall"), whether or not they fall within CN regions reporting deletions.
- OTV Caller is automatically run if any SNPs are assigned to the Conversion Type "OTV". Some of the resulting genotype calls may be "OTV", indicating a detected Off Target Variant.

With CN-aware Genotyping

- CN states will be reported for the predefined regions specified by the library package
- Unless overridden by the user, the control samples used for per-plate signal correction for CN analysis are identified by high concordance between measured and reference genotypes for expected control samples, and should also pass sample QC
- SNPs in a CN=0 region will be called as "ZeroCN"
- SNPs in a CN=1 region will be called with haploid genotype codes such as "A", "B", "NoCall_1"
- SNPs in a CN=2 or higher region will continue to be called with diploid genotype codes such as "AA", "AB", "BB", "NoCall"
- Special SNPs (those on chromosomes Y, MT, and non-PAR X) will also be assigned hemizygous and ZeroCN genotype codes when expected copy number is less than two
- CN-aware genotyping can be applied to probesets between two measured CN regions. For example, if two adjacent regions predict the same CN state, genotyping probesets between those two regions can be assigned genotype call codes consistent with that CN state.

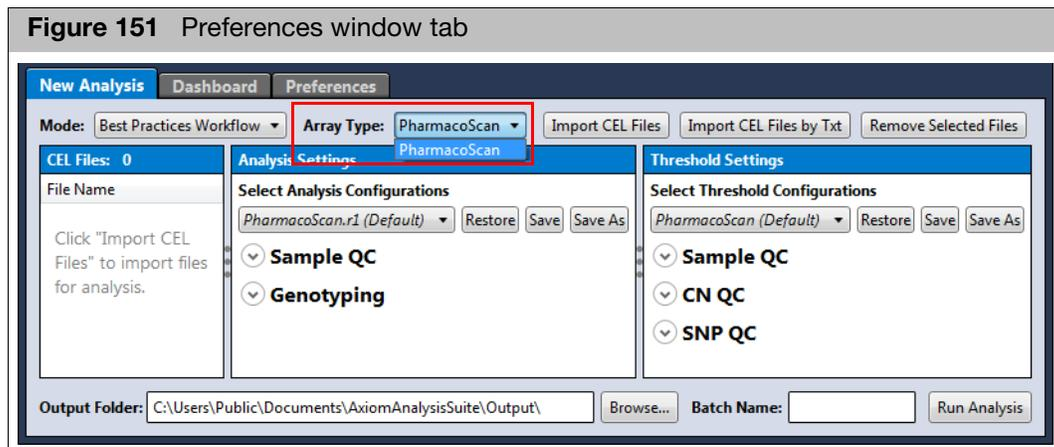
Note: Examples of library packages that support these copy number analyses include; CarrierScan for Predefined CN genotyping, and PharmacoScan for CN-aware genotyping.



Genotyping with CN Analysis

Setting Up a Genotyping with CN Analysis

Before setting up a genotyping with predefined CN analysis, click the New Analysis window tab, then click the Array Type drop-down menu to confirm an array such as PharmacoScan or CarrierScan is available, as shown in [Figure 151](#). If it is, continue to ["Selecting a Mode \(Workflow\)"](#).



Selecting a Mode (Workflow)

From the main Axiom Analysis Suite window tab, click the **Mode** drop-down.

1. Select **Best Practices Workflow**.
 Best Practices Workflow performs and combines the Sample QC Workflow and Genotyping Workflow. For more information, see ["Overview and Use of the Best Practices Workflow"](#) on page 137.

Importing CEL Files

1. Click **Import CEL Files**.
 The **Add CEL Files** window appears.
2. Navigate to your CEL file location.
3. Single-click on a CEL file or Ctrl click, Shift click, or press Ctrl A (to select multiple files).
4. Click **Open**.



The CEL Files pane populates and displays your selected CEL files. (Figure 152)

Figure 152 Cell File pane

CEL Files: 24	
File Name	
HG00127_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG00366_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG01770_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG02086_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG02301_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG02373_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG02589_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG02679_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03021_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03058_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03225_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03460_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03521_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03556_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03598_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
HG03643_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA18541_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA19118_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA19315_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA19318_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA19395_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA20289_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA20291_DB_10uLmPCRspike_Pharmacoscan_24_2016...	
NA20581_DB_10uLmPCRspike_Pharmacoscan_24_2016...	

After importing your CEL files, the Analysis and Threshold Setting panes auto-populate with default settings and additional user-configurable fields and settings specific to the library package, as shown in Figure 153.

For information on typical Axiom array analysis and threshold fields, see "Setting Up an Analysis" on page 19 and "Setting Up Threshold Settings" on page 28.

Figure 153 Example: Auto-populated Analysis and Threshold Setting panes

Analysis Settings

Select Analysis Configurations
 PharmacoScan_24F.r6 (Default) Restore Save Save As

Sample QC

GT Analysis File:
 PharmacoScan_24F_SNPSpecificPriors_Step1.r6.appt-genotype-axiom.AxiomGT1.appt2

Prior Model File:
 PharmacoScan_24F.r6.AxiomGT1

SNP List File:
 PharmacoScan_24F.r6.step1

Gender File (optional):

Hints/Inbred File (optional):
 Inbred Hints

Control Reference Calls File:
 PharmacoScan_24F.r6.signatureSNPs.refs.cn_controls

Genotyping

CN Control CEL List File (optional):

CN Analysis File:
 PharmacoScan_24F.r6.appt-copynumber-axiom-ssa.AxiomCN1.appt2

CN Reference Model File:
 PharmacoScan_24F.r6

CN Bins File:
 PharmacoScan_24F.r6.cnbins

GT Analysis File:
 PharmacoScan_24F_SNPSpecificPriors_Step2.r6.appt-genotype-axiom.AxiomGT1.appt2

Prior Model File:
 PharmacoScan_24F.r6.AxiomGT1

Multi-allele Background Prior Model File:
 PharmacoScan_24F.r6.AxiomGT1.mmb

Multi-allele Pairwise Prior Model File:
 PharmacoScan_24F.r6.AxiomGT1.mmp

Multi-allele Prior Model File:
 PharmacoScan_24F.r6.AxiomGT1.mm

SNP List File (optional):

Threshold Settings

Select Threshold Configurations
 PharmacoScan (Default) Restore Save Save As

Sample QC

Name	Settings
DQC	≥ 0.88
QC call_rate	≥ 98
Percent of passing samples	≥ 95
Average call rate for passing samples	≥ 98.5
Control comparisons	≥ 30
Control concordance	≥ 93

CN QC

Name	Settings
MAPD	≤ 0.35
Waviness SD	≤ 0.1

SNP QC

Name	Settings
species-type	Diploid
cr-cutoff	≥ 95
fld-cutoff	≥ 3.6
het-so-cutoff	≥ -0.2
het-so-XChr-cutoff	≥ -0.2
het-so-otv-cutoff	≥ -0.3
hom-ro-1-cutoff	≥ -0.5
hom-ro-2-cutoff	≥ -1.5
hom-ro-3-cutoff	≥ -1.5
hom-ro	true
hom-het	true
num-minor-allele-cutoff	≥ 2

Available for
CN-aware
genotyping
analysis variation

Available for
CN-aware
genotyping and
Predefined CN
analysis variations



Analysis Settings with Copy Number Options

X IMPORTANT! Only experienced users should modify default analysis settings.

Sample QC

- **GT Analysis File:** Parameters file for the genotyping step that calculates QC Call Rate.
- **Prior Model File:** Defines prior knowledge of SNP cluster locations. This file has the same format as a posteriors file, which is generated by the genotyping step. This means that you can "train" on a custom data set, and use the updated knowledge of cluster locations as a "seed" to possibly improve future genotyping batches. This file must contain two row entries for the GENERIC and GENERIC:1 probesets (if there are any probesets to be genotyped that are not listed in this file).
- **SNP List File:** A file of probeset IDs to genotype. For Sample QC it defines the probesets used to calculate QC Call Rate.
- **Gender File:** A file specifying the desired gender of every sample. If supplied, software will use values in this file instead of the computed gender. Gender impacts genotyping of chromosome X and Y SNPs.
- **Hints/Inbred File:** If a hints file, a file of expected genotype calls. This is used to influence the predicted cluster locations, which influences the final calls. In the Sample QC section, it influences the QC Call Rate.
- **Control Reference Calls File:** For CN-aware genotyping only, a recommended file containing the expected signature SNP calls of the CN control samples and used by Sample QC to identify the control samples among the supplied CEL files.

Genotyping

- **CN Control CEL List File:** This optional file identifies the CEL files that are to be used for per-plate tuning of CN signals. The file has the same format as file [Batch Name]\CNData\CNcontrolSamples.pass.txt, which is generated by a Sample QC step. The following table describes how the Workflow and this input option interact to select the controls used for CN analysis.

Workflow	CN Control CEL List File Input	Chosen Controls for CN Analysis
Best Practices	None	Passing controls identified by Sample QC step
Genotyping	None	No controls, therefore no plate signal tuning is done.
Best Practices	User-supplied	User-supplied controls
Genotyping	User-supplied	User-supplied controls

- **CN Analysis File:** Parameters file for the CN Analysis step.
- **CN Reference Model File:** Reference information for CN Analysis step.



- **CN Bins File:** Specifies for each predefined CN Region the callable CN states, and the MedianLog2Ratio boundaries defining each state.
- **GT Analysis File:** Parameters file for the final genotyping step.
- **Prior Model File:** Defines prior knowledge of bi-allelic SNP cluster locations. This file has the same format as a posteriors file, which is generated by the genotyping step. This means that you can “train” on a custom data set, and use the updated knowledge of cluster locations as a “seed” to possibly improve future genotyping batches. This file must contain two row entries for the GENERIC and GENERIC:1 probesets (if there are any probesets to be genotyped that are not listed in this file).
- If the library package supports multi-allele SNPs, then there are three additional model files for the three stages of multi-allele genotyping: Multi-allele Background Prior Model File, Multi-allele Pairwise Prior Model File, and Multi-allele Prior Model File.
- **SNP List File:** A file of probeset IDs to genotype. For Genotyping it defines the probesets for which genotypes will be reported.
- **Gender File:** A file specifying the known gender of every sample. If supplied, software will use values in this file instead of the computed gender. Gender impacts genotyping of chromosome X and Y SNPs.
- **Hints/Inbred File:** If a hints file, a file of expected genotype calls. This is used to influence the predicted cluster locations, which influences the final calls.
- **Posterior File Name:** The desired file output of a genotyping analysis, specifying identified SNP probeset cluster locations, variance, and relative weight. One use of this file is as a prior model file for future genotype analyses.
- **ps2snp File:** If multiple probeset designs exist on the array for a given SNP (for example, one forward and one reverse strand design), then the ps2snp file is used by the SNP classification step to identify the best performing probeset for the SNP, using the priority-order setting in the SNP QC section in the New Analysis tab. This text file has two tab delimited columns with the headers probeset_id and snpid (**snpid = affy_snp_id**).
- **Genotype Frequency File:** If the library package supports a check for unexpectedly high call frequency for specific genotypes, this optional file specifies the maximum expected frequency for reviewed genotypes.



Threshold Configurations with CN Options

✗ IMPORTANT! Only experienced users should modify default threshold settings.

Sample QC

Control Comparisons (CN-aware genotyping only): For identifying control samples, this is the minimum number of SignatureSNP probesets compared to a reference.

Control Concordance (CN-aware genotyping only): For identifying control samples, this is the minimum percent concordance of SignatureSNP calls to a reference.

CN QC

MAPD: Median Absolute Pairwise Difference of log₂ratio signals of adjacent copy number (CN) probesets must be below this value to make CN calls.

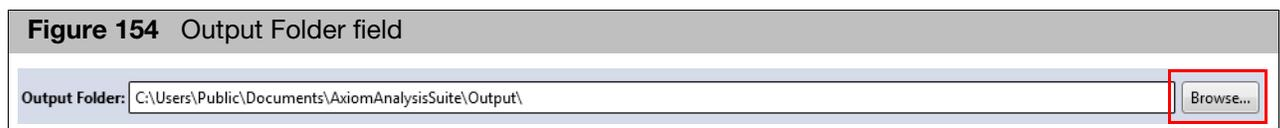
Waviness SD: Waviness Standard Deviation of log₂ ratio signals of copy number (CN) probesets must be below this value to make CN calls.

SNP QC

Refer to [Table 11 on page 149](#) for SNP QC Threshold name definitions.

Assigning an Output Folder Path

1. Click the **Output Folder** path's **Browse** button. ([Figure 154](#))



An Explorer window appears.

2. Navigate to the recommended path **C:\Users\Public\Documents\AxiomAnalysisSuite\Output**, then click **Select Folder**. Your selected output folder path is now displayed.

Assigning a Batch Name

The batch file is produced while your analysis is running and includes all the necessary files needed to view your analysis in the Axiom Analysis Suite Viewer.

1. Enter a name in the **Batch Name** field. ([Figure 155](#))



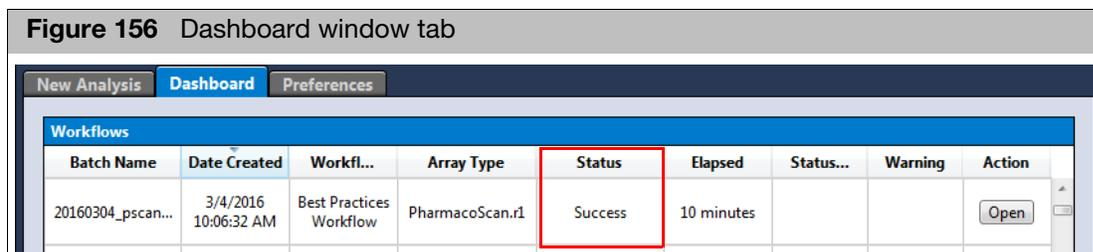
Note: A folder (with the same name as your entered batch name) is auto-generated during the analysis process. This folder includes all the necessary files needed to view your analysis results in the Viewer.



Running your Genotyping with CN Analysis

1. Click .

The Dashboard window tab appears. (Figure 156)



After **Success** is displayed in the Status column (Figure 156), click .

The button will not appear if a processing error occurs. If a message appears stating that no samples passed QC, click the **Open Selected Result(s)** button to review the analysis result that may need troubleshooting.

The Axiom Analysis Suite Viewer opens in a new window (Figure 157) and displays your completed analysis results.

Note: To make comparisons between your completed data analyses easier, open additional Viewer windows. To do this, click the **Dashboard** window tab, then click on the **Open** button again.

Viewing your Genotyping and CN Results

After processing arrays that support predefined CN or CN-aware genotyping, the following is added to the Axiom Analysis Suite Viewer:

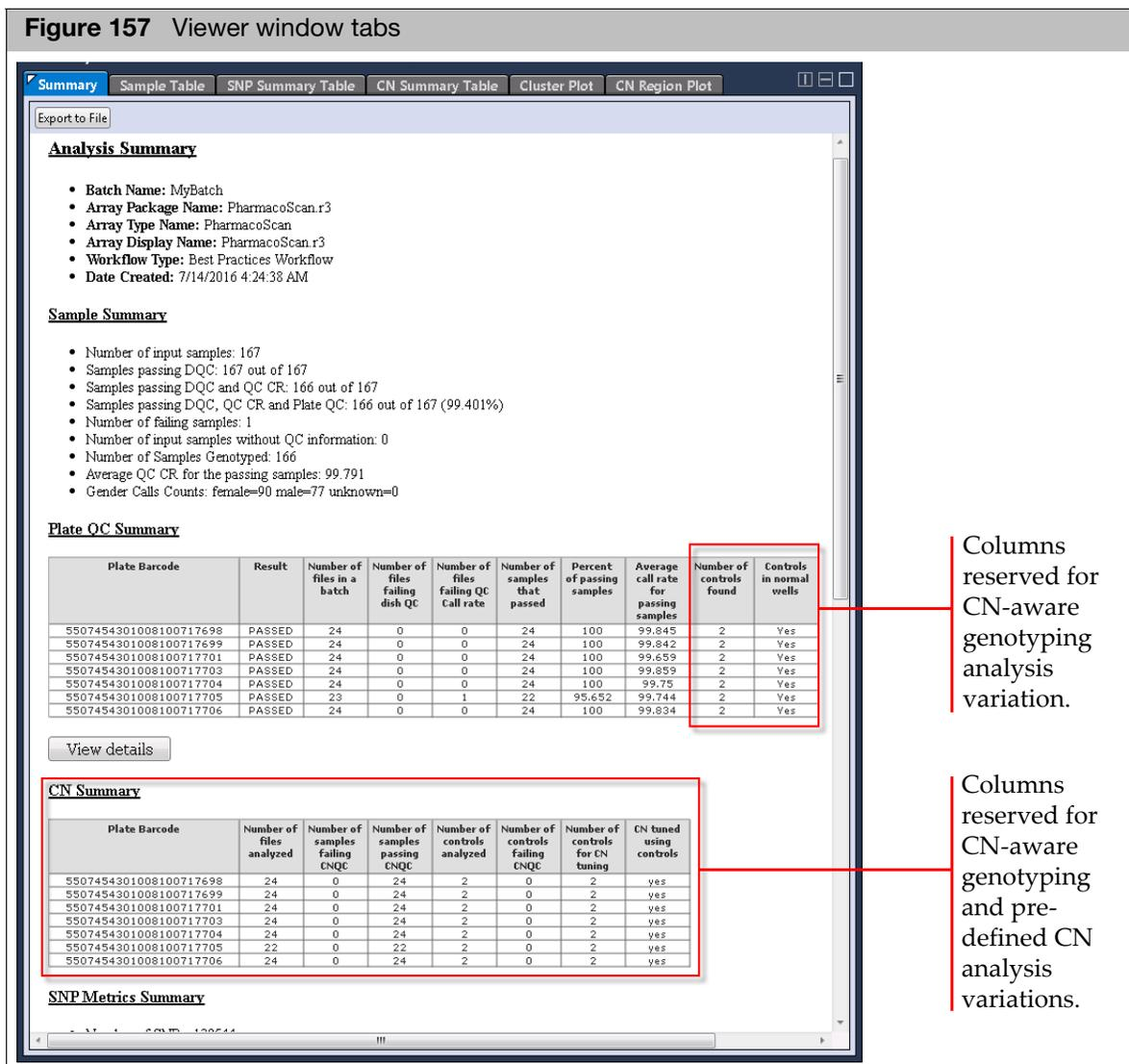
- New entries in the Summary Report, Sample Table, and SNP Summary Table.
- For CN-aware genotyping, the Cluster Plot also displays hemizygous and ZeroCN clusters for SNPs in CN regions.
- CN Summary Table, with Export CN Data option.
- A CN Region Plot window tab.
- Exported VCF formats can now include CN and SNP calls.



Summary Report

After successfully running a Sample QC or Best Practices workflow, the Summary report's **Plate QC Summary** section features two additional fields for the CN-aware genotyping analysis variation, as shown in [Figure 157](#).

- **Number of controls found** - This metric is a count of samples identified as controls based on high signature SNP concordance to expected calls in the Control Reference Calls library file.
- **Controls in normal wells** - The Control Reference Calls library file lists the expected plate wells for the control samples. This metric will report "Yes" if all the detected control samples are in the expected plate wells. It is OK to put the control samples in any wells you choose.



After successfully running a Genotyping or Best Practices workflow, the Summary report's **CN Summary** section features a table of information for each plate, as shown in [Figure 157](#). Refer to [Table 8](#) for descriptions of each CN Summary Report column.



Table 8 CN Summary Report

CN Summary Report	Description
Number of files analyzed	Count of CEL files supplied to the Genotyping step (which includes CN analysis).
Number of samples failing CNQC	Count of CEL files that fail copy number quality control checks, which is the result of a comparison of measured MAPD and waviness SD metrics against thresholds. Samples that fail CN QC report a CN state of NoCall. For the CN-aware genotyping analysis variation, SNPs from samples that fail CN QC are called without the benefit of CN-aware genotyping.
Number of samples passing CNQC	Count of CEL files that pass copy number quality control checks.
Number of controls analyzed	Count of CEL files supplied to the Genotyping step that are identified as controls.
Number of controls failing CNQC	<p>Count of control CEL files that fail copy number quality control checks, which is the result of a comparison of measured MAPD and waviness SD metrics against thresholds. Controls that fail CN QC report a CN state of NoCall. For the CN-aware genotyping analysis variation, SNPs from samples that fail CN QC are called without the benefit of CN-aware genotyping.</p> <p>Note: For the CN-aware genotyping analysis variation, Non-PAR X, Y, and mitochondrial SNPs use gender and other information to report non-diploid calls, regardless of CN QC status.</p>
Number of controls for CN tuning	Count of control CEL files that pass copy number quality control checks.
CN tuned using controls	If Number of controls for CN tuning > 0, then this value is "Yes". This means that CN signals are adjusted based on the signal measured in the control samples from the same plate. If the controls are appropriate for the samples being processed, this adjustment can improve CN calling accuracy.



Sample Table

After successfully running a CN analysis, additional columns appear within the Sample Table, as shown in [Figure 158](#).

Figure 158 Sample Table after a CN Analysis run

Sample Filename	Pass/Fail	DQC	QC call_rate	QC call_rate	QC het_rate	QC comput...	affymetrix-plate-barcode	affy... plate-peg...	Average call rate for passing samples	Percent of passing samples	MAPD	WavinessSD	CN passes MAPD	CN passes WavinessSD	CN passes QC	Control chosen	Control used	Number of controls for CN tuning	CN tuned using controls	
HG00185_JS...	Pass	0.978	99.935	99.85	29	27.415	male	5507454301008100717698	A05	99.844	100	0.264	0.043	yes	yes	yes	no	0	2	yes
HG00313_JS...	Pass	0.964	99.835	99.737	29.701	27.229	female	5507454301008100717698	C07	99.844	100	0.27	0.037	yes	yes	yes	no	0	2	yes
HG00404_JS...	Pass	0.959	99.825	99.812	25.298	23.805	female	5507454301008100717698	B05	99.844	100	0.291	0.033	yes	yes	yes	no	0	2	yes
HG00583_JS...	Pass	0.96	99.865	99.842	24.827	24.312	male	5507454301008100717698	D09	99.844	100	0.295	0.036	yes	yes	yes	no	0	2	yes
HG01085_JS...	Pass	0.969	99.87	99.805	29.316	28.885	male	5507454301008100717698	B09	99.844	100	0.258	0.041	yes	yes	yes	no	0	2	yes
HG01198_JS...	Pass	0.97	99.905	99.834	28.399	26.837	female	5507454301008100717698	E07	99.844	100	0.249	0.035	yes	yes	yes	no	0	2	yes
HG01519_JS...	Pass	0.974	99.915	99.843	29.396	27.048	female	5507454301008100717698	E05	99.844	100	0.308	0.043	yes	yes	yes	no	0	2	yes
HG02072_JS...	Pass	0.963	99.87	99.833	25.313	23.44	female	5507454301008100717698	B07	99.844	100	0.257	0.038	yes	yes	yes	no	0	2	yes
HG02086_JS...	Pass	0.956	99.87	99.77	25.649	23.964	female	5507454301008100717698	A09	99.844	100	0.265	0.035	yes	yes	yes	no	0	2	yes
HG02133_JS...	Pass	0.969	99.89	99.835	25.133	23.838	female	5507454301008100717698	E09	99.844	100	0.236	0.042	yes	yes	yes	no	0	2	yes
HG02589_JS...	Pass	0.956	99.78	99.751	24.336	23.467	female	5507454301008100717698	D07	99.844	100	0.27	0.03	yes	yes	yes	no	0	2	yes
HG02953_JS...	Pass	0.968	99.855	99.798	24.141	22.824	male	5507454301008100717698	F05	99.844	100	0.278	0.03	yes	yes	yes	no	0	2	yes
HG03460_JS...	Pass	0.958	99.82	99.725	24.056	22.894	male	5507454301008100717698	C05	99.844	100	0.326	0.036	yes	yes	yes	no	0	2	yes
NA18541_JS...	Pass	0.97	99.875	99.804	25.188	24.244	female	5507454301008100717698	D05	99.844	100	0.27	0.036	yes	yes	yes	no	0	2	yes
NA18642_T05...	Pass	0.969	99.835	99.782	25.393	24.289	female	5507454301008100717698	G05	99.844	100	0.254	0.036	yes	yes	yes	no	0	2	yes
NA19095_T04...	Pass	0.969	99.704	99.65	24.281	23.46	female	5507454301008100717698	H05	99.844	100	0.264	0.033	yes	yes	yes	no	0	2	yes
NA19114_JS...	Pass	0.968	99.9	99.763	23.805	23.175	female	5507454301008100717698	F07	99.844	100	0.241	0.03	yes	yes	yes	no	0	2	yes
NA19315_Ctrl...	Pass	0.962	99.915	99.84	24.281	23.732	female	5507454301008100717698	G09	99.844	100	0.254	0.024	yes	yes	yes	yes	1	2	yes
NA19318_Ctrl...	Pass	0.964	99.684	99.646	24.662	23.757	male	5507454301008100717698	H09	99.844	100	0.295	0.038	yes	yes	yes	yes	1	2	yes

Refer to [Table 9](#) for descriptions of each added Sample Table column.

Table 9 Added Sample Table columns after running a CN Analysis

Sample Table	Description
MAPD	Median Absolute Pairwise Difference of log2ratio signals of adjacent copy number (CN) probesets.
WavinessSD	Waviness Standard Deviation of log2ratio signals of adjacent copy number (CN) probesets
CN passes MAPD	"Yes" if the sample's MAPD value is not greater than the MAPD threshold used by CN QC.
CN passes WavinessSD	"Yes" if the sample's WavinessSD value is not greater than the WavinessSD threshold used by CN QC
CN passes QC	"Yes" if the sample passes both MAPD and WavinessSD threshold tests. If "no", the sample reports a CN state of NoCall. For the CN-aware genotyping analysis variation, SNPs from samples that fail CN QC are called without the benefit of CN-aware genotyping. Note: For the CN-aware genotyping analysis variation, Non-PAR X, Y, and mitochondrial SNPs use gender and other information to report non-diploid calls, regardless of CN QC status.
Control chosen	"Yes" if the control was listed in the CN Control CEL List file supplied for CN analysis.
Control used	"1" if the control was used to adjust CN signals of all samples in the same plate. The value will be 1 if Control chosen = 1 and CN passes QC = Yes.



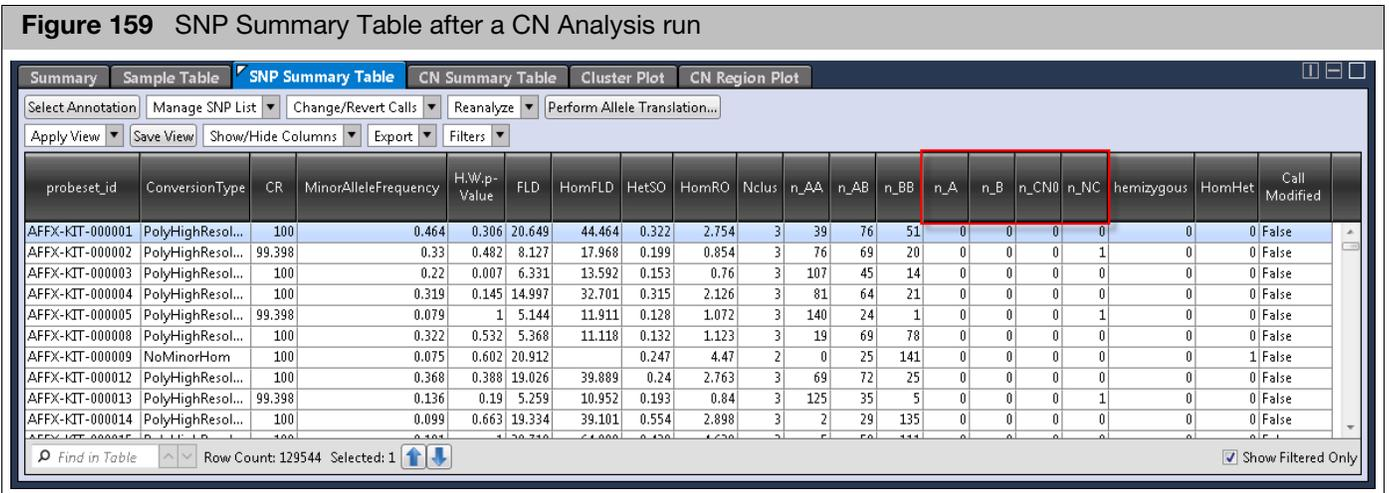
Table 9 Added Sample Table columns after running a CN Analysis

Sample Table	Description
Number of controls for CN tuning	A count of control samples used to adjust the CN signals for the given sample. This value will be the same for all samples on the same plate.
CN tuned using controls	If Number of controls for CN tuning > 0, then this value is "Yes". This means that CN signals are adjusted based on the signal measured in the control samples from the same plate. If the controls are appropriate for the samples being processed, this adjustment can improve CN calling accuracy.

SNP Summary Table

After successfully running a CN-aware genotyping analysis, additional columns appear within the SNP Summary Table, as shown in [Figure 159](#).

Figure 159 SNP Summary Table after a CN Analysis run



Refer to [Table 10](#) for descriptions of each added SNP Summary Table column.

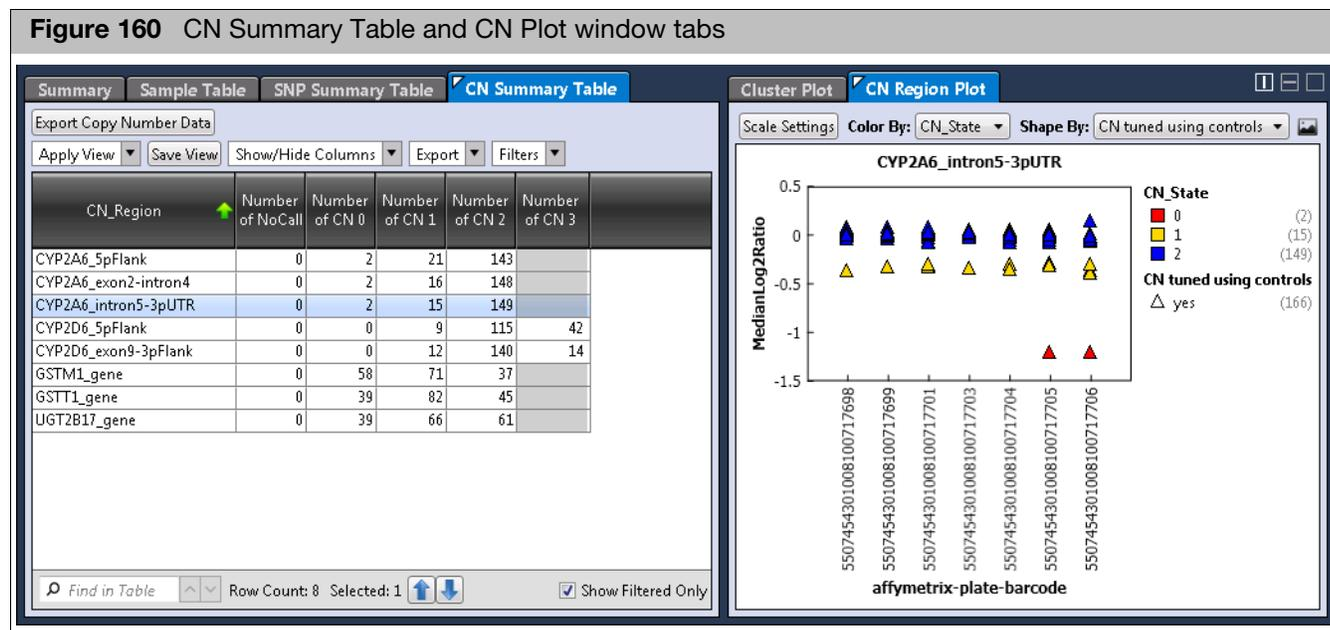
Table 10 Added SNP Summary Table columns after CN-aware genotyping is performed

SNP Summary Table	Description
n_A	Count of hemizygous "A" calls.
n_B	Count of hemizygous "B" calls.
n_CN0	Count of ZeroCN calls.
CopyNumIssue (hidden by default)	Copy Number Issue is "1" if the probeset is assigned to the "Other" ConversionType because the haploid or ZeroCN clusters are in an unusual location compared to each other or to the diploid clusters.



CN Summary Table and CN Region Plot

The Genotyping and Best Practices workflows perform predefined region copy number analysis (if the array content and library package support it). If this is the case, a **CN Summary Table** and **CN Region Plot** window tab appear in the Axiom Analysis Suite Viewer, as shown in [Figure 160](#).



CN Summary Table (Overview)

- The CN Summary Table gives a count of samples called with a given copy number state (CN_State) for a given copy number region (CN_Region).
- CN_States that cannot be reported are displayed as empty with a gray background. If the CN_State can be reported but no samples in the batch report that CN State, then the count is 0. As long as the default CN Bins library file is used during analysis setup, the only samples that will report a CN_State of NoCall are those that fail Copy Number QC.
- Selecting a specific CN_Region row in the table updates the associated CN Region Plot.
- The button **Export Copy Number Data** copies several CN reports to another location specified by the user. Note that CN_States for each CN_Region for each sample are also available in a VCF format export of genotyping data, available from the SNP Summary Table tab.



CN Region Plot (Overview)

- The CN Region Plot displays the MedianLog2Ratio of each genotyped sample for the copy number region selected in the CN Summary Table, grouped by affymetrix-plate-barcode.
- Selecting a sample or samples in the CN Region Plot also selects the same samples in the Sample Table and the Cluster Plot, and vice versa.
- Genes with a CN of 2 typically have a MedianLog2Ratio centered at 0, which indicates that the signals of probesets used for CN estimation are similar to the expected signals for a CN=2 sample.
- CN States are called by comparing the MedianLog2Ratio to the thresholds defined in the CN Bins library file selected during analysis setup. CN States cannot be edited in the Viewer.
- Sometimes between-plate assay variation can shift the observed MedianLog2Ratio values away from 0 for CN=2 samples. Fortunately, if the assay includes appropriate control samples on the assay plate, then the copy number signals can be tuned using the control samples. If CN plate correction was done, then the default "Shape By" metric "CN tuned using controls" identifies which samples had their CN measurements adjusted using the plate controls.

Overview and Use of the Best Practices Workflow

The recommended genotyping method is to use information from a batch of samples to improve the calling of individual samples. To this end, the cluster locations in signal space for each probeset (see the Cluster Plot) adapt in a Bayesian fashion to the supplied data. Prior knowledge of cluster locations influences the final calls. The more samples that are supplied in a batch, the more the final cluster locations will be influenced by the supplied data.

The dynamic nature of this genotyping algorithm means that if a given sample's CEL file is genotyped in a group of 24 CEL files, or in a group of 2400 CEL files, you can expect that some SNPs may experience a changed call (call <-> NoCall, or sometimes call 1 <-> call 2). This is more likely to happen if the SNP's data quality is not great (close clusters), if there are strong between-plate effects in cluster positions, or for very low Minor Allele Frequency SNPs where good information on rare genotype cluster positions may not be available before genotyping. Calling accuracy of both rare and common SNPs is improved even for small batch sizes if enough prior knowledge of cluster locations is available.

The batch nature of the genotyping means that if there are some samples of poor data quality in a group of good quality samples, then sometimes the poor quality samples will harm the calling accuracy of the good samples. For this reason, Axiom Analysis Suite's "Best Practices Workflow" can be thought of as a two-step process:

Step 1 Sample QC: Identify and exclude the poor quality sample CEL files.

Step 2 Genotyping: Genotype only the sample CEL files that pass Step 1.

- **Step 1:** The Sample QC Workflow performs the following steps:
 - Genotype a small set of SNPs used to uniquely identify each sample ("Signature SNPs"). This method uses static calling boundaries, so a given sample's calls are not influenced by other samples.
 - **DishQC:** Calculate DishQC metric and exclude CEL files with too small a DishQC from next step. For this method, a given sample's metrics are not influenced by other samples.
 - **QC Call Rate:** Initial cluster genotyping on remaining samples for selected QC SNPs, to identify and remove any additional CEL files with low QC call rates. For this method, batch information from other samples influences a given samples QC call rates. Genotypes are not stored.
 - **Plate QC:** Compare the average QC Call Rate of passing samples within a plate against the threshold "Average call rate for passing samples". If the metric is below this threshold, all the samples on the plate will fail Plate QC, and will not be genotyped.
 - **Identify Copy Number plate controls:**
 - For CN-aware genotyping arrays like PharmacoScan, auto-identify reagent control samples by comparing measured SignatureSNP genotypes from all CEL files against reference genotypes. A control is identified if [number of Signature SNPs with a call is \geq "Control comparisons"] AND [Concordance % of signature SNP calls compared with reference genotypes is \geq "Control concordance"]. Control samples that pass all sample QC checks will be used in the subsequent Genotyping workflow, unless overridden by the user.
 - For predefined CN region arrays like CarrierScan, select as CN controls all female samples passing sample QC.

- **Step 2:** Genotyping can be run by itself or as part of the Best Practices Workflow. When run as part of the Best Practices Workflow, only CEL files passing Sample QC are genotyped. When run by itself, the user has full control of which samples to genotype. Genotyping does the following steps:
 - For arrays supporting predefined region CN analysis or CN-aware genotyping, collate normalized signals for CN probesets of interest and perform predefined-region CN analysis. This step also calculates CN-specific QC metrics MAPD and wavinessSD. Samples that fail CN QC have their CN states reset to NoCall. If control samples are supplied (by user or by Sample QC step), then plate-based signal correction is performed. Control samples that fail CN QC are not used for plate-based signal correction.

X IMPORTANT! If you plan to run only the Genotyping workflow, and if the CN Control CEL List File is an Analysis Settings input option, this file should be supplied. If it is not, then plate signal correction is not performed and CN results may suffer.

The file **[Batch Name]\CNData\CNcontrolSamples.pass.txt** (from the Sample QC step) is supplied for CN analysis during a Best Practices Workflow.

- For CN-aware genotyping arrays like PharmacoScan, CN results are supplied as an input to the next genotyping step.
- Perform genotyping, using either default or user-supplied input options. Genotypes are saved.
- Generate summary statistics on each SNP.



About Allele Translation

Overview

Human genome sequence variation, which includes both single nucleotide polymorphisms (SNPs) as well as more complex structural variation in the form of insertions, duplications and deletions, underlies each individual's response to drugs. Products like PharmacoScan and DMET Plus are designed to enable comprehensive and accurate genotyping of specific polymorphisms involved in drug-metabolizing enzymes and transporters. Axiom Analysis Suite enables conversion of genotype calls to clinically-recognized star nomenclature via Allele Translation. This section explains the organization of the translation reports to help you interpret the translation data. Key concepts such as phase ambiguity and the impact of missing data on haplotype-based allele calling are described.

Gene Table Layout for Haplotyping

To appreciate how haplotyping operates, it is essential to describe the organization of the gene tables (Figure 161 and Figure 162) in the translation library file (*.translation).

Figure 161 Biological annotations [Example: Gene table data for markers in CYP1A1]

CYP1A1	Reference Link	Probe Set ID	Switch Design Strand to Report	dbSNP RS ID	Defining	cDNA		Change	Common Name
						Nucleotide Position	Genome Position		
CYP1A1	PMID: 11295847	AM_10774	Y	rs56313657	*6	993G>T	Ch15:75013804	M331I	CYP1A1*6_1635G>T(M331I)
CYP1A1	PMID: 15618738	AM_10771	Y	rs72547510	*7	1275_1276ir	Ch15:75013093	E426Fra	CYP1A1*7_2345insT
CYP1A1	PMID: 15618738	AM_10770	Y	rs72547509	*8	1343T>A	Ch15:75013026	I448N	CYP1A1*8_2413T>A(I448N)
CYP1A1	PMID: 8895751	AM_10769	Y	rs1799814	*4	1382C>A	Ch15:75012987	T461N	CYP1A1*4_2452C>A(T461N)
CYP1A1	PMID: 9070254	AM_10768	Y	rs1048943	*2C	1384A>G	Ch15:75012985	I462V	CYP1A1*2C_2454A>G(I462V)
CYP1A1	PMID: 11295847	AM_10766	N	rs41279188	*5	1390C>A	Ch15:75012979	R464S	CYP1A1*5or*9_2460C>A>T(R464SorC)
CYP1A1	PMID: 11295847	AM_10766	N	rs41279188	*9	1390C>T	Ch15:75012979	R464C	CYP1A1*5or*9_2460C>A>T(R464SorC)
CYP1A1	PMID: 15618738	AM_10765	Y	rs56240201	*10	1429C>T	Ch15:75012940	R477W	CYP1A1*10_2499C>T(R477W)
CYP1A1	PMID: 15618738	AM_10762	N	rs1800031	*3	*595T>C	Ch15:75012235	3'UTR	CYP1A1*3_3204T>C(3'UTR)
CYP1A1	PMID: 11295847	AM_10778	N	rs4646422	G45D	134G>A	Ch15:75015305	G45D	CYP1A1_134G>A(G45D)
CYP1A1	PMID: 11295847	AM_10776	N	rs34260157	R279W	835C>T	Ch15:75014049	R279W	CYP1A1_1390C>T(R279W)
CYP1A1	PMID: 9353182	AM_10775	Y	rs4987133	I286T	857T>C	Ch15:75014027	I286T	CYP1A1_1412T>C(I286T)
CYP1A1	PMID: 15618738	AM_10772	Y	rs2856833	F381L	1143C>A	Ch15:75013563	F381L	CYP1A1_1876C>A(F381L)
CYP1A1	GBID: NM_00049	AM_10767	Y	rs2278970	A463G	1388C>G	Ch15:75012981	A463G	CYP1A1_2458C>G(A463G)

Figure 162 Haplotype descriptions [Example gene table data for markers in CYP1A1]

Common Name	Haplotype	Reference Variant	*1	*2C	*3	*4	*5	*6	*7	*8	*9	*10	G45C	R279	I286T	F3
CYP1A1*6_1635G>T(M331I)	Y	G	T					T								
CYP1A1*7_2345insT	Y	-	T						T							
CYP1A1*8_2413T>A(I448N)	Y	T	A							A						
CYP1A1*4_2452C>A(T461N)	Y	C	A			A										
CYP1A1*2C_2454A>G(I462V)	Y	A	G		G											
CYP1A1*5or*9_2460C>A>T(R464SorC)	Y	C	A				A									
CYP1A1*5or*9_2460C>A>T(R464SorC)	Y	C	T							T						
CYP1A1*10_2493C>T(R477W)	Y	C	T								T					
CYP1A1*3_3204T>C(3'UTR)	Y	T	C		C											
CYP1A1_134G>A(G45D)	N	G	A										A			
CYP1A1_1390C>T(R279W)	N	C	T											T		
CYP1A1_1412T>C(I286T)	N	T	C												C	
CYP1A1_1876C>A(F381L)	N	C	A													A
CYP1A1_2458C>G(A463G)	N	C	G													

Biological annotations: The first set of columns in the table are annotations for the markers in each of the translated genes. Haplotype descriptions: The columns beyond the Common Name field contain information used for interpretation and translation of the gene file.

Following the columns enumerating the Reference and Variant alleles, haplotypes and markers in the gene are listed. In this example, CYP1A1 has 10 haplotypes described and they are named in the column headers. In addition, there are five additional markers for rare variants that can also be identified in this gene. Notice that the first haplotype described in this table is CYP1A1*1A, and that all markers except the last five are haplotyping markers (see the Haplotype field). Markers are characterized as "non-haplotyping" if their state in every reportable haplotype is not known. For example, a variant has been identified in this gene resulting in a non-synonymous change in the protein (CYP1A1_134G>A(G45D)), but the haplotype background of that variant is not available in the literature references used (for example, the Karolinska reference database; <http://www.cypalleles.ki.se/>).

Figure 161 illustrates the way that haplotypes are called. Notice that the haplotype names appear in the header of each gene table (for CYP1A1, when all markers are tested, this is *1, *2C, *3, *4, etc). Only differences from the reference haplotype appear in the haplotype columns, and the change for the altered base is indicated. For example, CYP1A1*2C contains one difference relative to CYP1A1*1, specifically a mutation 5'-prime to the gene, defined by the probe "CYP1A1*2C_2454A>G(I462V)".

The Defining field in the translation file lists the effect that a variant allele of this marker has. For example, because a variant at the final probe in the table, "CYP1A1_2458C>G(A463G)", results in a structural change in the protein, this marker is flagged with the name of the amino acid change that results: alanine at position 463 of the protein is a glycine in this variant (A463G). Although not shown in this gene, if a marker contains a variant allele in multiple haplotypes, then the marker does not uniquely define a single haplotype. That marker would then have an "N" (No) in the Defining field.

Haplotype field names with a # prefix, and rows with a # prefix in the Probe Set ID field are "commented out", and not used for translation.

Biological Annotations in Translation Reports

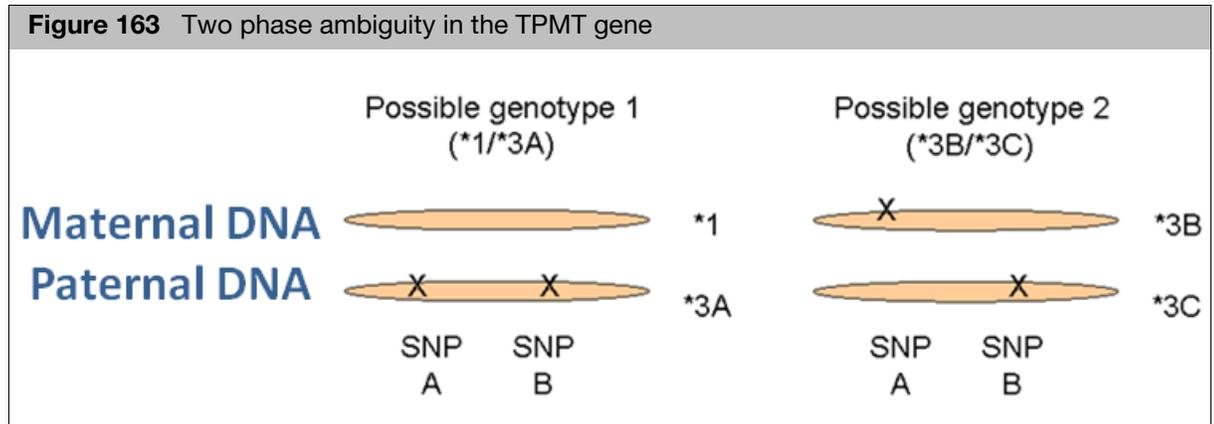
The primary function of the translation reporting is to summarize genotypes into commonly recognized variant names. In the case of the CYP450 core gene set, this translates to the Star-nomenclature followed by standardized nomenclature committee direction. Similar names are used by other steering committees such as the two Phase II enzyme genes, N-acetylase genes (NAT1 and NAT2) or the UGT-transferase gene families. Wherever possible, we have attempted to use a standard naming convention for the markers. To facilitate interpretation of the genotyping results, the translation reports provide:

- Reference publications, sequences or dbSNP identifiers for following previously published information about the variant site
- Precise genomic location in a recent genome build for identifying confirmatory genotyping assays
- Notation of protein changes that may result from the mutations in the panel. This field may also indicate whether the variant allele is strategically positioned in the promoter region or causes changes in splice junction sequences in the gene.
- Description of the initial star-allele which the variant was identified. Generally, this corresponds to the Summary flag entry
- Alternative alleles at each marker and whether the defined allele is the Reference base or Variant (corresponding to the altered gene form)

Along with the identified genotypes of the sample, this information provides biological evidence supporting haplotype calls.

Impact of Phase Ambiguity in Haplotyping

It is not unusual for individuals to be heterozygous at more than one defining marker (compound heterozygote genotype) in a gene. When this happens, multiple haplotype pairs may be consistent with resulting profiles, as shown in [Figure 163](#).



In this example, it is clear that the child could have inherited both variant alleles from one parent (in that case their diplotype would be *1/*3A), or could have inherited one variant allele from each of the two parents (in that case the diplotype would be *3B/*3C). Although the *3A haplotype is less common than either *3B or *3C, the translation reports list both potential haplotype pairs in the output reports. One reason for this is that the phenotypes may differ between the two alternative genetic configurations. In this case, three of the four alleles of TPMT have reduced activity: *3A, *3B and *3C, whereas the reference allele, *1 is a normally functioning allele. Phase ambiguity is relatively common in genes with common polymorphisms. [Figure 164 on page 143](#) lists the multiple possible calls due to phase ambiguity that were observed in six HapMap populations, and how often they occurred.

Figure 164 Example: Observed phase ambiguities in DMET Plus in a data set of six HapMap populations with 597 individuals and no children.

Gene	Call 1	Call 2	Call 3	Call Rate
CYP1A2	*1A/*1L	*1C/*1F		22.6%
CYP2B6	*1/*7	*5/*6		2.0%
CYP2D6	*2/*64	*10/*17		0.3%
FMO2	*2A/*3	*1/*2C		0.5%
NAT1	*4/*11	*11C/*30		1.8%
NAT2	*4/*5E	*5/*6		9.2%
NAT2	*4/*6J	*6/*7		3.7%
NAT2	*4/*14D	*6/*14		2.0%
NAT2	*4/*14F	*5/*14		1.3%
NAT2	*4/*7D	*7/*14		0.2%
SLC22A2	*1/*3D	*3A/*6		7.4%
SLC22A2	*1/*2B	*2A/*3A		6.2%
SLC22A2	*1/*3E	*2A/*3D	*2B/*6	5.5%
SLC22A2	*2B/*3D	*3A/*3E		1.3%
SLCO1B1	*1b/*17	*15/*21		5.9%
SLCO1B1	*1a/*14	*1b/*4		3.2%
SLCO1B1	*1a/*15	*1b/*5		3.2%
SLCO1B1	*1a/*17	*5/*21		2.5%
TPMT	*1/*3A	*3B/*3C		1.3%
UGT1A1	*1/*28+60	*28/*60		1.5%
UGT1A1	*1/*27+28+60+93	*27/*28+60+93		0.5%
UGT1ACOMMON	*76+79/*1A	*76/*79		8.0%
UGT2B15	*1/*5	*2/*4		12.9%

It is worth pointing out that the predicted phenotypes of some of these alternative diplotype calls are identical, and in these cases the Phenotype report will then report a single phenotype. When phase ambiguity is encountered and the Phenotype report does not resolve the multiple calls to a single Phenotype Call, follow-up metabolic screening may be merited to differentiate the actual genetic configuration of the test samples.

Diplotype to Phenotype Translation

Allele translations include a Phenotype report if the required metabolizer library file has been selected. The Phenotype report further translates the reported diplotypes (star allele pairs) from a subset of genes in the Comprehensive report into one of several phenotypes (e.g. "Poor Metabolizer"). As the software reads the comprehensive.rpt file, it will try to match Known Call diplotype values for each gene of each sample to one row of the metabolizer library file table. If a match is found, the associated phenotype and allele activities are written to the phenotype.rpt. If a match is not found, a Phenotype Call of "unknown" is reported. More information on this software feature is available in the DMET™ Plus Allele Translation white paper.

X IMPORTANT! Users are responsible for reviewing the metabolizer library file for accuracy!

Phenotype Call and Gene Activity interpretations for a Known Call are supported by differing levels of evidence from in vivo and/or in vitro research studies. Refer to metabolizer library file for a list of references. The actual phenotype and gene activities may be dependent on the substrate and dose.

If you do not want to generate a phenotype report, leave the Metabolizer File option blank in the Perform Allele Translation dialog.

If you want to report phenotypes for only a subset of genes, there are two ways to accomplish this:

1. Import a custom SNP List into the workspace containing probesets from only the genes of interest. At the point when you normally perform allele translation, select the option to filter to just probesets in this list.

Or

2. Create a copy of the metabolizer file that only contains the genes of interest.

If you want to change what phenotypes are reported for a particular combination of diplotypes, or you would like to change what is written to the header of the Phenotype Translation report, you will need to create and use a custom version of the .metabolizer library file. Instructions for doing this follow.

Creating a Custom Metabolizer Library File

If you choose to create a custom metabolizer table, start with the default metabolizer file as a template, then save a copy of this file using a new name.

X IMPORTANT! Use caution if editing the metabolizer file with Microsoft Excel. For example, Excel inserts quotation marks around text containing commas, which may make the file unreadable by Axiom Analysis Suite.

Before using the file with Axiom Analysis Suite, open it in another text editor and remove any unexpected text such as quotation marks.

To be recognized by Axiom Analysis Suite, the file:

- Must have the file extension ***.metabolizer**, where the ***** indicates your custom text.
- Must be encoded in ANSI, not Unicode or other encoding.
- Must exist in the library folder used by Axiom Analysis Suite.
- Must be selected for use from the **Configuration -> Options** menu, in the Translations tab.
- Must be properly formatted. It is recommended that you use a file comparison utility to verify that the only changes between the original and modified files are expected changes.

If you want to add phenotype reporting (for genes not currently in the metabolizer library file) the gene names and star allele names you want to add must exist in the ***.translation** library file used to generate the ***_comprehensive.rpt** file.

Metabolizer Library File Format

The **.metabolizer** library file is a tab-delimited text file that can be edited in any text editor. This file consists of a header section followed by a single table. Any rows from the start of the file until the beginning of the main table are considered header rows, and must begin with a pound or hash sign (#). Header rows are optional. Header rows beginning with **#%Info=** is added to the header of the ***_phenotype.rpt** file, so you can put custom text into your reports. The first row that does not begin with # must use the names shown below, be tab separated, and contain only lowercase letters.

```
gene allele_1 allele_2 phenotype activity_1 activity_2
```

Additional field names can be added to the first table row, but they will not be used. After the first table row, all rows require a value for the following fields:

```
gene allele_1 allele_2 phenotype
```

The **activity_1** and **activity_2** fields can be left blank. See [Figure 165 on page 147](#) for field descriptions.

Figure 165 field descriptions for the Metabolizer library file

Metabolizer field	Description														
gene	The gene name as reported in the comprehensive report. These values can also be found in the *.translation library file needed to generate the comprehensive report.														
allele_1 allele_2	The haplotype name of an allele for a gene as reported in the comprehensive report, e.g. '*2'. A Known Call in the comprehensive report is usually a single pair of alleles, e.g. '*1/*2'. To have this call be matched to a specific row in the metabolizer table, only one of the rows is needed in the following table: <table border="1" data-bbox="727 583 1321 800"> <thead> <tr> <th>gene</th> <th>allele_1</th> <th>allele_2</th> <th>phenotype</th> </tr> </thead> <tbody> <tr> <td>CYP2D6</td> <td>*1</td> <td>*2</td> <td>NM</td> </tr> <tr> <td>CYP2D6</td> <td>*2</td> <td>*1</td> <td>NM</td> </tr> </tbody> </table> <p>Axiom Analysis Suite will report an error IF it detects duplicate rows (as in the above example) AND IF the duplicate rows report disagreeing phenotypes.</p>	gene	allele_1	allele_2	phenotype	CYP2D6	*1	*2	NM	CYP2D6	*2	*1	NM		
gene	allele_1	allele_2	phenotype												
CYP2D6	*1	*2	NM												
CYP2D6	*2	*1	NM												
phenotype	The value that should be reported for the associated 'gene allele_1/allele_2' call. The string should be short with no commas, quotes, or whitespace characters. Common phenotype names are: <table border="1" data-bbox="727 1024 1247 1520"> <thead> <tr> <th>phenotype</th> <th>definition</th> </tr> </thead> <tbody> <tr> <td>UM</td> <td>Ultra-rapid metabolizer</td> </tr> <tr> <td>RM</td> <td>Rapid metabolizer</td> </tr> <tr> <td>NM</td> <td>Normal metabolizer</td> </tr> <tr> <td>IM</td> <td>Intermediate metabolizer</td> </tr> <tr> <td>PM</td> <td>Poor metabolizer</td> </tr> <tr> <td>unknown</td> <td>Unknown metabolizer</td> </tr> </tbody> </table>	phenotype	definition	UM	Ultra-rapid metabolizer	RM	Rapid metabolizer	NM	Normal metabolizer	IM	Intermediate metabolizer	PM	Poor metabolizer	unknown	Unknown metabolizer
phenotype	definition														
UM	Ultra-rapid metabolizer														
RM	Rapid metabolizer														
NM	Normal metabolizer														
IM	Intermediate metabolizer														
PM	Poor metabolizer														
unknown	Unknown metabolizer														
activity_1 activity_2	The reported gene activity or function level for an allele, e.g. 'normal' or 'decreased'. Activity_1 is for allele_1, and activity_2 is for allele_2. The values in these fields are used to populate the Gene Activity field in the phenotype report, e.g. 'normal/decreased'. If you leave these fields empty, the phenotype report will display '/' for the Gene Activity.														
Optional fields	Axiom Analysis Suite will ignore additional fields in the metabolizer file. Additional fields may be used to annotate each row.														

Reference Databases Used in Translation Data Curation

The databases used to curate the allele translation gene tables include:

- PharmGKB - Stanford University Pharmacogenomics reference database
<http://www.pharmgkb.org>
- Karolinska cytochrome P450 gene standard nomenclature
<http://www.cypalleles.ki.se>
- Database of NAT genes (Democritus University of Thrace)
<http://nat.mbg.duth.gr>
- Database of UGT genes
<https://www.pharmacogenomics.pha.ulaval.ca/ugt-alleles-nomenclature>
- Drug interaction database (University of Indiana)
<http://medicine.iupui.edu/clinpharm/ddis>
- PubMed - On-line National Library of Medicine publication database
<http://www.ncbi.nlm.nih.gov/pubmed>



Threshold Names

Use the table of definitions below to help select thresholds (when setting up an analysis or regenerating SNP metrics).

Table 11 Threshold Names

Threshold Group	Threshold Name	Description
Sample QC	DQC	A sample's dish QC value must be greater than or equal to 'DQC' to pass sample QC.
Sample QC	QC call_rate	A sample's call rate value must be greater than or equal to 'QC call_rate' to pass genotyping QC.
Sample QC	Percent of passing samples	If a plate's percent of passing samples is smaller than this number, all samples on the plate will show a warning in the Sample Table.
Sample QC	Average call rate for passing samples	A plate's average QC call rate of passing samples must be greater than or equal to this number to pass plate QC.
Sample QC	Control comparisons	CN-aware genotyping parameter. For identifying control samples, this is the minimum number of SignatureSNP probesets compared to a reference.
Sample QC	Control concordance	CN-aware genotyping parameter. For identifying control samples, this is the minimum percent concordance of SignatureSNP calls to a reference.
CN QC	MAPD	For arrays that support copy number (CN) analysis, the Median Absolute Pairwise Difference of log ₂ ratio signals of adjacent copy number (CN) probesets must be less than this value to make CN calls.
CN QC	Waviness SD	For arrays that support copy number (CN) analysis, the Waviness Standard Deviation of log ₂ ratio signals of copy number (CN) probesets must be less than this value to make CN calls.
SNP QC	species-type	Species type for the array, which affects some SNP QC checks. Species types include: Diploid, Human, and Polyploid. Note: Selecting 'Human' will assign a probeset with two clusters that are both homozygous to a non-recommended ConversionType like 'Other'.
SNP QC	cr-cutoff	Minimum acceptable call rate.

**Table 11** Threshold Names

Threshold Group	Threshold Name	Description
SNP QC	fld-cutoff	For autosomal probesets, minimum acceptable FLD value for cluster separation.
SNP QC	het-so-cutoff	Minimum acceptable value for the correctness of the Size (Y position) offset of the heterozygous cluster.
SNP QC	het-so-XChr-cutoff	For probesets on the non-pseudoautosomal regions of chromosome X, the minimum acceptable value for the correctness of the Size (Y position) offset of the female heterozygous cluster.
SNP QC	het-so-otv-cutoff	Minimum acceptable value for the correctness of the Size (Y position) offset of the heterozygous cluster, possibly indicating a fourth cluster below the heterozygous cluster (OTV).
SNP QC	hom-ro-1-cutoff	Minimum acceptable value for the correctness of the Contrast (X position) of the homozygous clusters (Ratio Offset) when a probeset has 1 genotype cluster.
SNP QC	hom-ro-2-cutoff	Minimum acceptable value for the correctness of the Contrast (X position) of the homozygous clusters (Ratio Offset) when a probeset has 2 genotype clusters.
SNP QC	hom-ro-3-cutoff	Minimum acceptable value for the correctness of the Contrast (X position) of the homozygous clusters (Ratio Offset) when a probeset has 3 genotype clusters.
SNP QC	hom-ro	Flag indicating whether the metric HomRO is used in classification.
SNP QC	hom-het	If hom-het flag is FALSE, a probeset with two clusters that are homozygous and heterozygous is assigned a non-recommended ConversionType like 'Other'. Hom-het flag should be TRUE for human and diploid species for a large batch of samples, and should be FALSE for highly inbred samples such as polyploid species.
SNP QC	num-minor-allele-cutoff	Minimum minor allele count for categorizing a probeset as PolyHighResolution.
SNP QC	hom-ro-hap-1-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 1 haploid genotype cluster.
SNP QC	hom-ro-hap-1-XChr-cutoff	For non-PAR X probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 1 haploid genotype cluster.
SNP QC	hom-ro-hap-1-YChr-cutoff	For Y probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 1 haploid genotype cluster.

**Table 11** Threshold Names

Threshold Group	Threshold Name	Description
SNP QC	hom-ro-hap-1-MTChr-cutoff	For mitochondrial probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 1 haploid genotype cluster.
SNP QC	hom-ro-hap-2-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 2 haploid genotype clusters.
SNP QC	hom-ro-hap-2-XChr-cutoff	For non-PAR X probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 2 haploid genotype clusters.
SNP QC	hom-ro-hap-2-YChr-cutoff	For Y probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 2 haploid genotype clusters.
SNP QC	hom-ro-hap-2-MTChr-cutoff	For mitochondrial probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters from 0 (Ratio Offset) when a probeset has 2 haploid genotype clusters.
SNP QC	hom-hap-X-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Contrast (X position) of the haploid clusters relative to the homozygous clusters.
SNP QC	hom-hap-Y-lower-cutoff	For autosomal probesets, maximum acceptable value for the correctness of the Size (Y position) of the haploid clusters relative to homozygous clusters.
SNP QC	hom-hap-Y-upper-cutoff	For autosomal probesets, maximum acceptable value for the correctness of the Size (Y position) of the haploid clusters relative to homozygous clusters.
SNP QC	CN0-hap-X-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Contrast (X position) of the ZeroCN cluster relative to the haploid clusters.
SNP QC	CN0-hap-X-YChr-cutoff	For Y probesets, minimum acceptable value for the correctness of the Contrast (X position) of the ZeroCN cluster relative to the haploid clusters.
SNP QC	CN0-hap-Y-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Size (Y position) of the ZeroCN cluster relative to the haploid clusters.
SNP QC	CN0-hap-Y-YChr-cutoff	For Y probesets, minimum acceptable value for the correctness of the Size (Y position) of the ZeroCN cluster relative to the haploid clusters.
SNP QC	CN0-dip-X-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Contrast (X position) of the ZeroCN cluster relative to the diploid clusters.

**Table 11** Threshold Names

Threshold Group	Threshold Name	Description
SNP QC	CN0-dip-Y-cutoff	For autosomal probesets, minimum acceptable value for the correctness of the Size (Y position) of the ZeroCN cluster relative to the diploid clusters.
SNP QC	aaf-XChr-cut	For non-PAR X probesets, maximum acceptable difference in male versus female A-allele frequency. Any probeset that fails is categorized as Other.
SNP QC	fld-XChr-cut	For non-PAR X probesets, minimum acceptable FLD value for diploid cluster separation. Note: FLD is computed differently for non-PAR X probesets than for autosomal probesets.
SNP QC	homfld-XChr-cut	For non-PAR X probesets, minimum acceptable FLD value for male haploid cluster separation.
SNP QC	homfld-YChr-cut	For Y probesets, minimum acceptable FLD value for male haploid cluster separation.
SNP QC	sign-diff-hom-1-cutoff	For multi-allele probesets, minimum acceptable difference between mean log2 signal and background when there is 1 homozygous genotype cluster.
SNP QC	sign-diff-hom-2-cutoff	For multi-allele probesets, minimum acceptable difference between mean log2 signal and background when there are 2 homozygous genotype clusters.
SNP QC	min-mean-cp2-cutoff	For multi-allele probesets, minimum acceptable mean log2 allele signal for samples that appear to have at least 2 copies of that allele.
SNP QC	max-mean-cp2-cutoff	For multi-allele probesets, maximum acceptable mean log2 allele signal for samples that appear to have at least 2 copies of that allele.
SNP QC	min-genotype-freq-samples	Minimum count of samples to support the genotype frequency p-value calculation. This parameter is used if a genotype frequency file is supplied.
SNP QC	genotype-p-value-cutoff	Minimum acceptable value for the genotype frequency p-value calculation. Probesets not meeting this threshold may be categorized as 'UnexpectedGenotypeFreq'. This parameter is used if a genotype frequency file is supplied, and if the count of genotyped samples is at least min-genotype-freq-samples.
SNP QC	priority-order	Priority order of probeset conversion types when performing probeset selection.
SNP QC	recommended	Probesets having ConversionTypes in this recommended set will be included in the recommended probeset list.
SNP QC	y-restrict	Y-restrict is the maximum vertical distance 2 points in 2 clusters can be when matching up points for calculating the edge metric values BB_dis_x_adj and AA_dis_x_adj. Y-restrict is not used to set probeset ConversionType. It is only used during 'Regenerate SNP Metrics' when the 'Generate advanced metrics' option is selected.



Table 11 Threshold Names

Threshold Group	Threshold Name	Description
PS Supplemental	homfld-cut	Minimum acceptable FLD value for homozygous cluster separation.
PS Supplemental	bb-varx-cut	Maximum acceptable Contrast (X position) variance for BB cluster. Not used if bb-varx-z-cut-enabled is true.
PS Supplemental	bb-vary-cut	Maximum acceptable Size (Y position) variance for BB cluster. Not used if bb-vary-z-cut-enabled is true.
PS Supplemental	ab-varx-cut	Maximum acceptable Contrast (X position) variance for AB cluster. Not used if ab-varx-z-cut-enabled is true.
PS Supplemental	ab-vary-cut	Maximum acceptable Size (Y position) variance for AB cluster. Not used if ab-vary-z-cut-enabled is true.
PS Supplemental	aa-varx-cut	Maximum acceptable Contrast (X position) variance for AA cluster. Not used if aa-varx-z-cut-enabled is true.
PS Supplemental	aa-vary-cut	Maximum acceptable Size (Y position) variance for AA cluster. Not used if aa-vary-z-cut-enabled is true.
PS Supplemental	bb-varx-z-cut-enabled	If true, bb-varx-z-cut is used.
PS Supplemental	bb-vary-z-cut-enabled	If true, bb-vary-z-cut is used.
PS Supplemental	ab-varx-z-cut-enabled	If true, ab-varx-z-cut is used.
PS Supplemental	ab-vary-z-cut-enabled	If true, ab-vary-z-cut is used.
PS Supplemental	aa-varx-z-cut-enabled	If true, aa-varx-z-cut is used.
PS Supplemental	aa-vary-z-cut-enabled	If true, aa-vary-z-cut is used.
PS Supplemental	bb-varx-z-cut	Maximum acceptable Contrast (X position) Z-score for BB cluster. Not used if bb-varx-z-cut-enabled is false.
PS Supplemental	bb-vary-z-cut	Maximum acceptable Size (Y position) Z-score for BB cluster. Not used if bb-vary-z-cut-enabled is false.
PS Supplemental	ab-varx-z-cut	Maximum acceptable Contrast (X position) Z-score for AB cluster. Not used if ab-varx-z-cut-enabled is false.
PS Supplemental	ab-vary-z-cut	Maximum acceptable Size (Y position) Z-score for AB cluster. Not used if ab-vary-z-cut-enabled is false.
PS Supplemental	aa-varx-z-cut	Maximum acceptable Contrast (X position) Z-score for AA cluster. Not used if aa-varx-z-cut-enabled is false.
PS Supplemental	aa-vary-z-cut	Maximum acceptable Size (Y position) Z-score for AA cluster. Not used if aa-vary-z-cut-enabled is false.
PS Supplemental	clustermin	Minimum number of samples in a cluster for the new variance or Z-score value to be calculated.
PS Supplemental	n-minor-hom-cut	Minimum number of minor hom samples for PolyHighResolution classification. PolyHighResolution probesets where the number of minor hom samples is less than n-minor-hom-cut are categorized as nMinorHom.

**Table 11** Threshold Names

Threshold Group	Threshold Name	Description
PS Supplemental	hetv-maf-cut	Maximum acceptable heterozygosity parameter. PolyHighResolution probesets with het rate > MAF*hetv-maf-cut are categorized as UnexpectedHeterozygosity.
PS Supplemental	variance-class	List of categories that supplemental filters are used on.

Sample Table

Use the table of definitions below for the fields in the Sample Table.

Table 12 Sample Table

Selection	Definition
Sample Filename	CEL file name.
Pass/Fail	Sample quality control status. A sample is called 'Pass' by the Sample QC step if it meets the minimum thresholds for DQC, QC call_rate, and average call rate for passing samples.
DQC	DishQC measures the amount of overlap between two homozygous peaks created by non-polymorphic probes. DQC of 1 is no overlap, which is good. DQC of 0 is complete overlap, which is bad.
call_rate	Percentage of autosomal SNPs with a call other than NoCall. 'SpecialSNP' probesets are excluded, as they are mostly non-autosomal.
QC call_rate	Percentage of autosomal SNPs with a call other than NoCall (measured at the Sample QC step).
het_rate	Percentage of SNPs called AB (i.e. the heterozygosity) for autosomal SNPs.
QC het_rate	Percentage of SNPs called AB (i.e. the heterozygosity) for autosomal SNPs (measured at the Sample QC step).
computed_gender	Computed gender for the sample.
QC computed_gender	Computed gender for the sample (measured at the Sample QC step).
affymetrix-plate-barcode	Plate barcode number.
QC affymetrix-plate-barcode	Plate barcode number (measured at the Sample QC step).
affymetrix-plate-peg-wellposition	Well position of the plate's peg.
QC affymetrix-plate-peg-wellposition	Well position of the plate's peg (measured at the Sample QC step).
Average call rate for passing samples	Average QC Call Rate of passing samples within the plate to which this sample belongs.
Percent of passing samples	Percentage of samples passing sample QC within the plate to which this sample belongs.
cel_filepath	CEL file path.
cel_file_identifier	CEL file identifier.

**Table 12** Sample Table

Selection	Definition
affymetrix-array-id	Command Console array identifier.
total_call_rate	Call rate at the default or user-specified threshold for all SNPs.
QC total_call_rate	Call rate at the default or user-specified threshold for all SNPs (measured at the Sample QC step).
total_het_rate	Percentage of SNPs called AB (i.e., the heterozygosity) for all SNPs.
QC total_het_rate	Percentage of SNPs called AB (i.e., the heterozygosity) for all SNPs (measured at the Sample QC step).
hom_rate	Percentage of SNPs called AA or BB (i.e. the homozygosity) for autosomal SNPs.
QC hom_rate	Percentage of SNPs called AA or BB (i.e. the homozygosity) for autosomal SNPs (measured at the Sample QC step).
total_hom_rate	Percentage of SNPs called AA or BB (i.e. the homozygosity) for all SNPs.
QC total_hom_rate	Percentage of SNPs called AA or BB (i.e. the homozygosity) for all SNPs (measured at the Sample QC step).
cluster_distance_mean	Average distance to the cluster center for the called genotype.
QC cluster_distance_mean	Average distance to the cluster center for the called genotype (measured at the Sample QC step).
cluster_distance_stdev	Standard deviation of the distance to the cluster center for the called genotype.
QC cluster_distance_stdev	Standard deviation of the distance to the cluster center for the called genotype (measured at the Sample QC step).
allele_summarization_mean	Average of the allele signal estimates (log ₂ scale).
QC allele_summarization_mean	Average of the allele signal estimates (log ₂ scale) (measured at the Sample QC step).
allele_summarization_stdev	Standard deviation of the allele signal estimates (log ₂ scale).
QC allele_summarization_stdev	Standard deviation of the allele signal estimates (log ₂ scale) (measured at the Sample QC step).
allele_deviation_mean	Average of the absolute difference between the log ₂ allele signal estimate and its median across all arrays.
QC allele_deviation_mean	Average of the absolute difference between the log ₂ allele signal estimate and its median across all arrays (measured at the Sample QC step).
allele_deviation_stdev	Standard deviation of the absolute difference between the log ₂ allele signal estimate and its median across all arrays.
QC allele_deviation_stdev	Standard deviation of the absolute difference between the log ₂ allele signal estimate and its median across all arrays (measured at the Sample QC step).
allele_mad_residuals_mean	Average of the median absolute deviation (MAD) between observed probe intensities and probe intensities fitted by the model.

**Table 12** Sample Table

Selection	Definition
QC allele_mad_residuals_mean	Average of the median absolute deviation (MAD) between observed probe intensities and probe intensities fitted by the model (measured at the Sample QC step).
allele_mad_residuals_stdev	Standard deviation of the median absolute deviation (MAD) between observed probe intensities and probe intensities fitted by the model.
QC allele_mad_residuals_stdev	Standard deviation of the median absolute deviation (MAD) between observed probe intensities and probe intensities fitted by the model (measured at the Sample QC step).
A_signal_mean	Average of control A probe raw intensities in the AT channel.
T_signal_mean	Average of control T probe raw intensities in the AT channel.
G_signal_mean	Average of control G probe raw intensities in the GC channel.
C_signal_mean	Average of control C probe raw intensities in the GC channel.
AT_B	AT channel background, which is the average signal of the GC control probes in AT channel.
GC_B	GC channel background, which is the average signal of the AT control probes in GC channel.
AT_S	AT channel signal, which is the average signal of the AT control probes in AT channel.
GC_S	GC channel signal, which is the average signal of the GC control probes in GC channel.
AT_SBR	AT channel signal to background ratio, defined as AT_S/AT_B.
GC_SBR	GC channel signal to background ratio, defined as GC_S/GC_B.
AT_B_IQR	AT channel interquartile range (middle 50%) of background intensities, measured using GC control probes.
GC_B_IQR	GC channel interquartile range (middle 50%) of background intensities, measured using AT control probes.
AT_S_IQR	AT channel interquartile range (middle 50%) of signal intensities, measured using AT control probes.
GC_S_IQR	GC channel interquartile range (middle 50%) of signal intensities, measured using GC control probes.
CV_AT	AT channel median coefficient of variation of replicate probe signals for control probesets.
CV_GC	GC channel median coefficient of variation of replicate probe signals for control probesets.
AT_FLD	Fisher's Linear Discriminant between signal and background in the AT channel, defined as $[\text{median_of_AT_probe_intensities} - \text{median_of_GC_probe_intensities}]^2 / [0.5 * (\text{AT_S_IQR}^2 + \text{AT_B_IQR}^2)]$.



Table 12 Sample Table

Selection	Definition
GC_FLD	Fisher's Linear Discriminant between signal and background in the GC channel, defined as $[\text{median_of_GC_probe_intensities} - \text{median_of_AT_probe_intensities}]^2 / [0.5 * (\text{GC_S_IQR}^2 + \text{GC_B_IQR}^2)]$.
log_diff_qc	A cross channel QC metric, defined as $\text{mean}(\log(\text{AT_SBR})) / \text{std}(\log(\text{AT_SBR})) + \text{mean}(\log(\text{GC_SBR})) / \text{std}(\log(\text{GC_SBR}))$, where signal and background are calculated for control non-polymorphic probes after intensity normalization.
saturation_AT	Fraction of features in the AT channel with intensity greater than or equal to 3800. Features likely to saturate the scanner in the long exposure image will instead be measured in the short exposure image, and all feature signals are scaled using a high dynamic range exposure merging technique.
saturation_GC	Fraction of features in the GC channel with intensity greater than or equal to 3800. Features likely to saturate the scanner in the long exposure image will instead be measured in the short exposure image, and all feature signals are scaled using a high dynamic range exposure merging technique.
cn-probe-chrXY-ratio_gender_meanX	Average probe intensity (raw, untransformed) of X chromosome nonpolymorphic probes.
cn-probe-chrXY-ratio_gender_meanY	Average probe intensity (raw, untransformed) of Y chromosome nonpolymorphic probes.
cn-probe-chrXY-ratio_gender_ratio	Gender ratio $Y/X = \text{cn-probe-chrXY-ratio_gender_meanY} / \text{cn-probe-chrXY-ratio_gender_meanX}$.
cn-probe-chrXY-ratio_gender	Predicted gender, based on the value of cn-probe-chrXY-ratio_gender_ratio .
reagent_version	Reagent version used for processing the arrays, based on data intensity values. Note: You can only perform batch genotyping analysis on CEL files processed using the same reagent version.
reagent_discrimination_value	Value assigned to the reagent, and used to determine reagent_version .
user_color	User-supplied color to associate with the sample. For more information, see " Setting User Colors " on page 55.



Concordance Columns

Use the table of definitions below for the columns in the Concordance table.

Table 13 Concordance Columns

Column Name	Description
Base	The first sample in the comparison.
Reference	The second sample in the comparison.
#SNPs Called	Number of SNPs common to both sample and reference files with genotype calls.
#Concordant SNP's	Number of called SNPs that have the same genotype call.
%Concordance	Percentage of called SNPs that have the same genotype call.



Annotations and Columns

Use the table of definitions below to help select the annotations you can add to the SNP Summary Table.

Table 14 Annotations and Columns

Column Name	Description
Probe Set ID	The Affymetrix unique identifier for the set of probes used to detect a particular Single Nucleotide Polymorphism (SNP probe sets only).
Affx SNP ID	The Affymetrix unique identifier for the set of probes used to detect a particular Single Nucleotide Polymorphism (SNP). (SNP probe sets only, not available for Axiom™ Genome-Wide Human Array).
dbSNP RS ID	The dbSNP ID that corresponds to this probe set or SNP. The dbSNP at the National Center for Biotechnology Information (NCBI) attempts to maintain a unified and comprehensive view of known single nucleotide polymorphisms (SNPs), small scale insertions/deletions, polymorphic repetitive elements, and microsatellites from TSC and other sources. The dbSNP is updated periodically, and the dbSNP version used for mapping is given in the dbSNP version field. For more information, see: http://www.ncbi.nlm.nih.gov/SNP/ (SNP probe sets only).
Chromosome	The chromosome on which the SNP is located on the current Genome Version.
Chromosome Start	The nucleotide base start position where the SNP is found. The genomic coordinates given are in relation to the current genome version and may shift as subsequent genome builds are released.
Chromosome Stop	The nucleotide base stop position where the SNP is found. The genomic coordinates given are in relation to the current genome version and may shift as subsequent genome builds are released.
Strand	Genomic strand that the SNP resides on.
Cytoband	Cytoband location of the SNP derived from the SNP physical map and the chromosome band data provided by UCSC.
Strand Vs dbSNP	Indicates whether the SNP is on the same or reverse strand as compared to dbSNP (SNP probe sets only).
ChrX pseudo-autosomal region	SNPs on the X Chromosome which are mapped to the two pseudo-autosomal region have a value of 1 or 2 in this field. All other SNPs are indicated by 0. A value of "1" indicates that the marker maps to the PAR-1 region and a value of "2" indicates that the marker maps to the PAR-2 region. A value of "0" indicates that the marker does not map to either of the two PAR regions.
Probe Count	The total number of probes in the probe set.
Flank	The nucleotide sequence surrounding the SNP. This is a 33-mer sequence with 16 nucleotides on either end of the SNP position. The alleles at the SNP position are provided in the brackets (SNP probe sets only).

**Table 14** Annotations and Columns

Column Name	Description
Allele A, Allele B	<p>At array (or underlying database) design time, the following naming convention is used to assign allele nucleotide bases to the "Abstract" allele codes "A" and "B":</p> <ol style="list-style-type: none"> 1. SNPs are fixed on the forward strand of the design-time reference genome. 2. For AT or CG SNPs (SNP alleles are A/T or C/G), the alleles are named in alphabetical order (A and C are the "A" alleles, in these cases); 3. For non-AT and non-CG SNPs, allele A is A or T, allele B is C or G; 4. For indels, allele A is -, allele B is the insertion. 5. For multi-base alleles, the alleles are named in alphabetical order. (For [AGT/TTA], AGT would be "Allele A". For [GGT/TTA], GGT would be "Allele A".)
Ref Allele, Alt Allele	<p>The reference allele and alternative alleles are specified according to the current reference genome build. The value of Ref Allele could be "-", which indicates an insertion after the specified position. Otherwise it is the sequence of the allele of the marker which matches the current reference genome. The value of Alt Allele could be "-", which indicates that the variant is a deletion with respect to the current genome build. Otherwise it is the sequence(s) of the allele(s) of the marker which does not match the current reference genome. If neither allele of the marker matches the current genome build sequence, then the value of the Ref Allele is set to "." and the value of the Alt Allele(s) is set to <allele_1>/<allele_2>, where <allele_1> and <allele_2> are the alleles of the marker. If the current genome build position of the marker is unknown then the value of the Ref Allele and Alt Allele are set to "---" to denote missing information.</p>
Associated Gene	<p>SNPs were associated with human genes by comparing the genomic locations of the SNPs to genomic alignments of human mRNA sequences. In cases where the SNP is within a known gene, NetAffx reports the association. Additionally, for genes with exon or CDS annotations, NetAffx reports whether or not the SNP is in an exon, and in the coding region. If the SNP is not within a known gene, NetAffx reports the closest genes in the genomic sequence, and the distance and relationship of the SNP relative to the genes. A SNP is upstream of a gene if it is located closer to the 5' end of the gene and is downstream of a gene if it is located closer to the 3' end of the gene.</p>
Genetic Map	<p>Describes the genetic location of the SNP derived from three separate linkage maps (deCODE, Marshfield, or SLM). The physical distance between the markers is assumed to be linear with their genetic distance. The genetic location is computed using the linkage maps from the latest physical location of the SNP and the neighboring microsatellite markers (SNP probe sets only).</p>
Microsatellite	<p>Describes the nearest microsatellite markers (upstream, downstream and overlapping) for the SNP.</p>
Enzyme Fragment	<p>Lists the enzyme, the restriction fragment containing the SNP and the fragment length. The Whole Genome Assay protocol detects SNPs that are contained within the genomic restriction fragments to simplify the sequence background for genotyping arrays (not available for Axiom Genome-Wide Human Array).</p>
Copy Number Variation	<p>When available, a description of Copy Number Variation Region (CN) probe sets as described by the Database of Genomic Variants (not available for Axiom Genome-Wide Human Array).</p>

**Table 14** Annotations and Columns

Column Name	Description
SNP Interference	This column is for Copy Number probe sets. It indicates whether or not a known SNP overlaps a copy number probe (CN probe sets only, not available for Axiom Genome-Wide Human Array).
In Final List	This column annotates extended content for genotyping arrays. A value of "1" indicates that the marker is included in the final version of the library file and a value of "0" indicates that the marker is not included in the final version of the library file (SNP probe sets only, not available for Axiom Genome-Wide Human Array).
% GC	The fraction of bases that are G or C in a window of 250,000 bases to each side of the SNP or CN position. All positions that are nearer to the end than 250,001 are set to the value of the position at 250,001 from that end. Position and chromosome values for SNPs and CN probes were mapped to the position of bases in the FASTA files for the build of the genome used in this release of NetAffx, and these bases were then used for all calculations (not available for Axiom Genome-Wide Human Array).
Heterozygous Allele Frequencies	Describes the heterozygous frequency of the allele from Yoruba, Japanese, Han Chinese and CEPH studies using the Affymetrix genotyping arrays. (SNP probe sets only)
Allele Sample Size	Sample size used for Allele Frequency estimates (SNP probe sets only).
Allele Frequencies	Describes the major and minor frequency of the allele from Yoruba, Japanese, Han Chinese and CEPH studies using the Affymetrix genotyping arrays (SNP probe sets only).
Minor Allele	Indicates the Minor Allele of a SNP (SNP probe sets only).
Minor Allele Frequency	The Minor Allele Frequency of a SNP (SNP probe sets only).
OMIM ID	Furnishes OMIM and Morbid Map IDs and their respective gene titles. This database contains information from the Online Mendelian Inheritance in Man® (OMIM®) database, which has been obtained under a license from the Johns Hopkins University. This database/product does not represent the entire, unmodified OMIM® database, which is available in its entirety at www.ncbi.nlm.nih.gov/omim/ .
Ordered Alleles	A list of alleles alphabetically ordered by abstract allele code. For bi-allele probesets, the order is "Allele A // Allele B" For multi-allele probesets, the order is "Allele A // Allele B // Allele C // Allele D", with the number of alleles equal to Allele Count.
Allele Count	The number of measured alleles. For bi-allele probesets the count is 2. For multi-allele probesets, the count is 3 or more.



SNP Summary Table Definitions

Use the table of definitions below to help select your SNP Summary Table columns.

Note: Most SNP Summary Table metrics are calculated only for the subset of samples reported by the column 'gender_metrics'. Refer to the gender_metrics description for more information.

Note: Some of the metrics defined in the following table are not displayed by default in the SNP Summary Table. Select **Show/Hide Columns** or **Apply View > All Columns View** to access more metrics. Some metrics are reported only after selecting the option(s) **Generate advanced metrics** or **Run PS Supplemental**, when choosing to **Regenerate SNP Metrics**. Some metrics are reported only if multi-allele probesets are present, or if CN-aware genotyping was performed.

Note: A few metrics are calculated only for multi-allele probesets. Most metrics are calculated only for biallele probesets.

Table 15 SNP Summary Table Metrics

Column Name	Description
ConversionType	Probeset classification
CR	Call rate (CR) is the percentage of samples with a genotype call other than "No Call" for the SNP. Note: Call Rate on non-pseudoautosomal regions of chromosome X is reported only for female samples. Call Rate on chromosome Y is reported only for male samples.
MinorAlleleFrequency	<p>The allele frequency for the A allele is calculated as:</p> $P_A = \frac{(\# AA \text{ Calls} + 0.5 * AB \text{ Calls})}{\text{Total \# Calls}}$ <p>Where the Total # Calls does not include the No Calls. The B allele frequency is .</p> $P_B = 1 - P_A$ <p>The minor allele frequency is the Min(P_A, P_B).</p>



Table 15 SNP Summary Table Metrics

Column Name	Description
<p>H.W.p-Value</p>	<p>Hardy Weinberg p-value is a measure of the significance of the discrepancy between the observed ratio or heterozygote calls in a population and the ratio expected if the population was in Hardy Weinberg equilibrium.</p> <p>There are two statistical tests used for HWE. When AA, AB, and BB counts are all ≥ 10, a Chi-squared test is used. When one or more of the AA, AB, and BB counts are < 10, an Exact test is used. An Exact test means that the p-value is calculated exactly and not approximated from a population distribution.</p> $x^2 = \frac{(f^2_{aa} - f_a)^2}{f^2_{aa}} + \frac{(2f_{aa}f_{bb} - f_{ab})^2}{2f_{aa}f_{bb}} + \frac{(f^2_{bb} - f_b)^2}{f^2_{bb}}$ <p>Where:</p> $f_a = \frac{(\#AA \text{ Calls})}{\text{Total \# Calls}}$ $f_b = \frac{(\#BB \text{ Calls})}{\text{Total \# Calls}}$ $f_{aa} = \frac{(\#AA \text{ Calls} + 0.5 * \# AB \text{ Calls})}{\text{Total \# Calls}}$ $f_{bb} = \frac{(\#BB \text{ Calls} + 0.5 * \# AB \text{ Calls})}{\text{Total \# Calls}}$ $f_{ab} = \frac{(\#AB \text{ Calls})}{\text{Total \# Calls}}$ $PHW = CDF(x^2)$ <p>Where CDF is the Cumulative Distributive Function for the chi-squared distribution.</p> <p>The Exact test used is the one implemented in R package "HardyWeinberg" for more information see:</p> <p>[2] Haldane, J., 1954. An exact test for randomness of mating. J. Genet. 52 631-635.</p> <p>[3] Levene, H., 1949. On a matching problem arising in genetics. Ann. Math. Stat. 20 91-94.</p>
<p>H.W.statistic</p>	<p>H.W.statistic is 1 if H.W.p-Value is calculated using an exact test. H.W.statistic is 0 if the chi-squared test is used. Refer to H.W.p-Value definition (above) for more information.</p>
<p>minGenotypeFreqPval</p>	<p>The minimum probability that any of the clusters present for a probeset have more observations than the supplied expected values. If the ConversionType is 'UnexpectedGenotypeFrequency' then minGenotypeFreqPval is smaller than the SNP QC parameter genotype-p-value-cutoff. The metric minGenotypeFreqPval is available for supported arrays if a genotype frequency file is supplied for SNP QC, and if the genotyped sample batch size is at least as large as the SNP QC parameter 'min-genotype-freq-samples'.</p>

**Table 15** SNP Summary Table Metrics

Column Name	Description
minGenotypeFreqCluster	The genotype cluster for which minGenotypeFreqPval is reported.
FLD	Fisher's Linear Discriminant (FLD) is a measure of the cluster quality of a probeset. High-quality probeset clusters have well-separated centers, and the clusters are narrow. FLD is measured from the SNP posteriors produced during genotyping. FLD is undefined if either the heterozygous or one of the homozygous clusters is empty.
HomFLD	HomFLD is a version of FLD computed for the homozygous genotype clusters. HomFLD is undefined for probesets without two homozygous clusters.
HomFLD_hap	HomFLD_hap is a version of HomFLD computed for the haploid genotype clusters, but only for samples specified by gender_metrics. HomFLD_hap is undefined for probesets without two haploid clusters.
HetSO	Heterozygous Strength Offset measures how far the heterozygous cluster center sits above the homozygous cluster centers in the Size dimension (Y position). Low HetSO values are produced either by misclustering events or by the inclusion of samples that contain variations from the reference genome. Most well-clustered diploid SNPs have positive HetSO values.
HomRO	Homozygote Ratio Offset is the distance to zero in the Contrast dimension (X position) from the center of the homozygous cluster that is closest to zero. If there is only one homozygous cluster, HomRO is the distance from that cluster center to zero on the Contrast dimension.
HomRO_hap	HomRO_hap is a version of HomRO computed for haploid clusters, but only for samples specified by gender_metrics. HomRO_hap will report -10 otherwise.
nMinorAllele	The count of minor alleles, which is one for each heterozygous call, one for each minor haploid call, and two for each minor homozygous call.
Nclus	The number of genotype clusters.
n_AA	The number of AA calls.
n_AB	The number of AB calls.
n_BB	The number of BB calls.
n_A	The number of A calls (haploid).
n_B	The number of B calls (haploid).
n_CN0	The number of ZeroCN calls (zero copy number).
n_NC	The number of NoCall calls, including NoCall_1 (haploid).
AA.meanX	Average Contrast (X position) for AA cluster.
AA.meanY	Average Size (Y position) for AA cluster.
AB.meanX	Average Contrast (X position) for AB cluster.
AB.meanY	Average Size (Y position) for AB cluster.
BB.meanX	Average Contrast (X position) for BB cluster.

**Table 15** SNP Summary Table Metrics

Column Name	Description
BB.meanY	Average Size (Y position) for BB cluster.
hemizygous	Hemizygous flag is 1 if the probeset measures chromosome Y or mitochondrial DNA, indicating that diploid genotypes are not possible. Otherwise the flag is 0.
BB_dis_x_adj	Smallest Contrast (X position) distance between any BB cluster sample and any neighbor cluster sample, adjusted to only comparing the sample pairs that have a smaller Size (Y position) distance than y_restrict parameter.
AA_dis_x_adj	Smallest Contrast (X position) distance between any AA cluster sample and any neighbor cluster sample, adjusted to only comparing the sample pairs that have a smaller Size (Y position) distance than y_restrict parameter.
freq_diff	A-allele frequency difference between male and female samples.
HomHet	HomHet flag is 1 if, when two diploid genotype clusters are present, one cluster is homozygous and the other is heterozygous. Otherwise the flag is 0.
BB.varX	Contrast (X position) variance for BB cluster.
BB.varY	Size (Y position) variance for BB cluster.
AB.varX	Contrast (X position) variance for AB cluster.
AB.varY	Size (Y position) variance for AB cluster.
AA.varX	Contrast (X position) variance for AA cluster.
AA.varY	Size (Y position) variance for AA cluster.
BB.varX.Z	Contrast (X position) variance Z-score for BB cluster.
BB.varY.Z	Size (Y position) variance Z-score for BB cluster.
AB.varX.Z	Contrast (X position) variance Z-score for AB cluster.
AB.varY.Z	Size (Y position) variance Z-score for AB cluster.
AA.var.X.Z	Contrast (X position) variance Z-score for AA cluster.
AA.var.Y.Z	Size (Y position) variance Z-score for AA cluster.
CopyNumIssue	CopyNumIssue flag is 1 if the probeset is categorized as 'Other' because of unusual locations of the haploid or ZeroCN clusters. Otherwise the flag is 0.
BestProbeset	BestProbeset flag is available when multiple probesets are mapped to the same SNP (Affy-SNP-ID) by a ps2snp file. A probeset is selected based on the priority order of the conversion types. BestProbeset flag is 1 when it is the best or only probeset for a SNP. Otherwise the flag is 0.
BestandRecommended	BestandRecommended flag is 1 if BestProbeset is 1 and the ConversionType belongs to the Recommended set of conversion types. Otherwise the flag is 0.

**Table 15** SNP Summary Table Metrics

Column Name	Description
gender_metrics	Not all probeset metrics are reported using all samples. Number of clusters (Nclus) is calculated on all samples. See the Call Rate (CR) definition for samples used. The remaining metrics are calculated using the sample set described in gender_metrics. If gender_metrics is 'all', all samples are used. If gender_metrics is 'female', only female samples are used. If gender_metrics is 'male', only male samples are used. If gender_metrics is 'diploid', then haploid and ZeroCN samples are excluded before calculation.
MinMean_cp_2	For multi-allele probesets, the minimum mean log2 allele signal for samples that appear to have at least two copies of that allele. If this value is not between the thresholds for the SNP QC parameters 'min-mean-cp2-cutoff' and 'max-mean-cp2-cutoff', then the probeset fails this QC test.
minSigBgndDiffHom	For multi-allele probesets, minimum difference between mean log2 signal and background for the homozygous genotype clusters. If this value is less than the threshold for the appropriate SNP QC parameter 'sign-diff-hom-1-cutoff' or 'sign-diff-hom-2-cutoff', then the probeset fails this QC test.
count_ma_A, count_ma_B, count_ma_C, ...	For multi-allele probesets, the count of the associated allele. For example, if there are only two samples with the calls AA and AC, then count_ma_A is 3 and count_ma_C is 1.
Call Modified	Call Modified flag is True if any calls for this probeset are changed since the batch results were first created. If no calls are changed the flag is False.

For support visit [thermofisher.com/support](https://www.thermofisher.com/support) or email techsupport@lifetech.com
thermofisher.com

18 August 2017

ThermoFisher
SCIENTIFIC