



User Manual

Somatic Mutation Viewer

For Research Use Only. Not for use in diagnostic procedures.

Trademarks

Affymetrix®, GeneChip®, NetAffx®, Command Console®, Powered by Affymetrix™, GeneChip-compatible™, Genotyping Console™, DMET™, GeneTitan®, Axiom®, CytoScan®, and GeneAtlas® are trademarks or registered trademarks of Affymetrix, Inc. All other trademarks are the property of their respective owners.

All other trademarks are the property of their respective owners.

This database/product 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 <http://www.omim.org/>

Limited License Notice

Limited License. Subject to the Affymetrix terms and conditions that govern your use of Affymetrix products, Affymetrix grants you a non-exclusive, non-transferable, non-sublicensable license to use this Affymetrix product only in accordance with the manual and written instructions provided by Affymetrix. You understand and agree that except as expressly set forth in the Affymetrix terms and conditions, that no right or license to any patent or other intellectual property owned or licensable by Affymetrix is conveyed or implied by this Affymetrix product. In particular, no right or license is conveyed or implied to use this Affymetrix product in combination with a product not provided, licensed or specifically recommended by Affymetrix for such use.

Patents

Software products may be covered by one or more of the following patents: U.S. Patent Nos. 5,733,729; 5,795,716; 5,974,164; 6,066,454; 6,090,555; 6,185,561; 6,188,783; 6,223,127; 6,228,593; 6,229,911; 6,242,180; 6,308,170; 6,361,937; 6,420,108; 6,484,183; 6,505,125; 6,510,391; 6,532,462; 6,546,340; 6,687,692; 6,607,887; 7,062,092 and other U.S. or foreign patents.

Copyright

© 2013 Affymetrix, Inc. All rights reserved.

Contents

Chapter 1	Somatic Mutation Viewer	2
	Introduction	2
	System Requirements	2
	Installing Somatic Mutation Viewer	2
	Starting Somatic Mutation Viewer	2
<hr/>		
Chapter 2	Setting Up the Viewer	4
	Loading OSCHP Files	4
	Assigning a Reference Model File	5
	Assigning an Annotation File	6
<hr/>		
Chapter 3	Using the Viewer	7
	OSCHP Files Window	7
	OSCHP File Window Columns	7
	Adding and Removing OSCHP Files	8
	Sorting OSCHP File Window Columns	8
	Sample View Tab Window	9
	Using the Sample View Tab Window	9
	MutScores (Scores) and MutCalls (Calls) - Overview	11
	Marker Information Window	12
	Editing Thresholds in the Marker Information Window	13
	Marker View Tab Window	15
	Editing Thresholds in the Marker View Tab Window	15
	Saving your New Thresholds	16
	Viewing Tools	17
	Resizing Window Panes	17
	Taskbar Options	17
	Box Whisker Plot Graphic	18
	Saving your Edited OSCHP Files	18

Somatic Mutation Viewer

Introduction

This document provides the general guidelines on how to use the Somatic Mutation Viewer application.

Somatic Mutation Viewer is a tool to visualize somatic mutation MutScores and MutCalls across a set of supplied OncoScan FFPE Assay OSCHP files.

In addition to visualizing the MutScores, the tool also enables you to edit the thresholds used to make the MutCalls and save its updated calls to OSCHP files.

You can view the data grouped by marker or by sample, and you can select individual data points for more information. Summary statistics are provided by sample or by marker. Sort these tables by the field of interest to quickly find rows of interest.

System Requirements

Operating System
Windows® 7 Professional (64-bit) with Service Pack 1 installed

Installing Somatic Mutation Viewer

To install Somatic Mutation Viewer:

1. Go to **www.affymetrix.com** and navigate to the Somatic Mutation Viewer location.
2. Locate and download the zipped Somatic Mutation Viewer software package.
3. Unzip the file, then double-click **SomaticMutationApp.exe** to install it.
4. Follow the directions provided by the installer.

Starting Somatic Mutation Viewer

To start Somatic Mutation Viewer:

1. Locate the Somatic Mutation Viewer Shortcut  on your system's Desktop, then double-click on it.

The Viewer opens. (Figure 1.1 on page 3)

Figure 1.1 Somatic Mutation Viewer- Main window

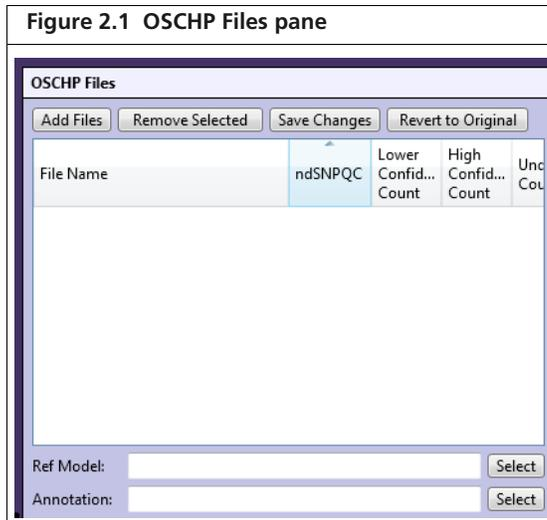
The Somatic Mutation Viewer main window displays the following components:

- Header:** affymetrix logo and Somatic Mutation Viewer title.
- Tabs:** Sample View (selected) and Marker View.
- Plot Area:** A large empty plot area with seven vertical lines representing markers. The values for these markers are: -1.94, -0.97, 0.00, 0.97, 1.94, 2.91, and 3.87.
- Bottom Left Controls:** Min: -2, Max: 5, Reset Scale, Copy to Clipboard.
- OSCHP Files Panel:**
 - Buttons: Add Files, Remove Selected, Save Changes, Revert to Original.
 - Table with columns: File Name, ndSNPQC, Lower Confid... Count, High Confid... Count, Uncou...
 - Ref Model: [Select]
 - Annotation: [Select]
- Marker Information Panel:**
 - Table with columns: Probeset Name, Low Thresh..., High Thresh..., Lower Confid... Count, High Confid... Count.
 - Buttons: Load Thresholds, Create Threshold File, Reset Thresholds.

Setting Up the Viewer

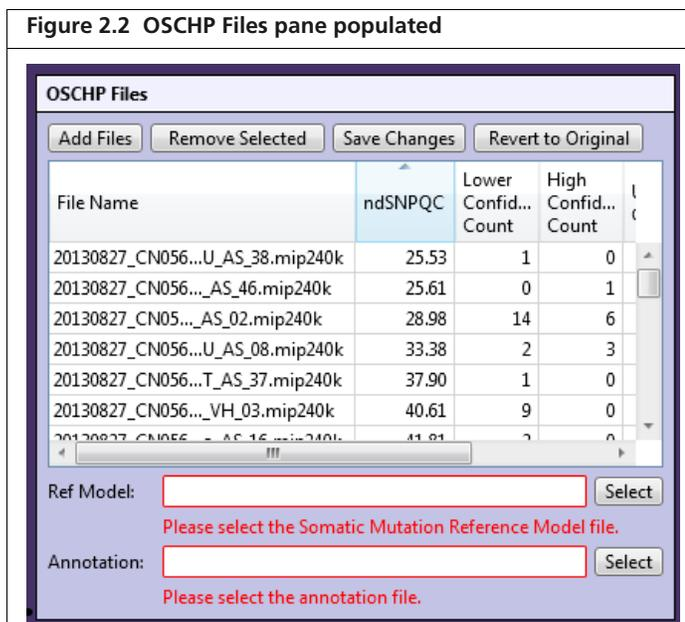
IMPORTANT: The Somatic Mutation Viewer requires OncoScan array data (OSCHP files).

Loading OSCHP Files

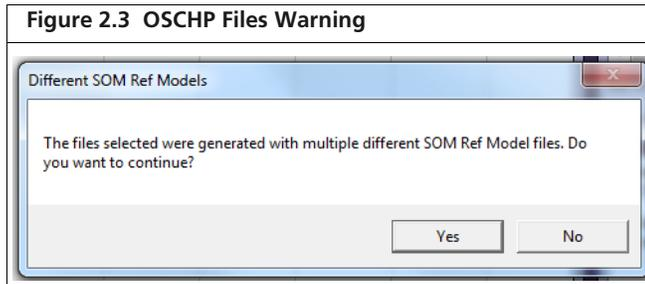


To Load OSCHP Files:

1. Click **Add Files**.
A File window appears.
2. Navigate to your OSCHP folder, then single click, Ctrl click, Shift click, or Ctrl-A (to select multiple or all OSCHP files).
3. Click **Open**.
The OSCHP File Name pane is now populated. (Figure 2.2)



A warning may appear if the selected OSCHP files were generated with different SOM Ref Model File. (Figure 2.3)



Do one of the following:

- Click **Yes** to continue.
- Click **No** to remove the current list of OSCHP files.

Assigning a Reference Model File



NOTE: The reference model file supplies the information that generates the Marker View's ox Box Whisker Plot graphic. Make sure to select the same reference model file you used to generate the OSCHP data.

To select a Reference Model file:

1. Click **Select**.

A File window appears.

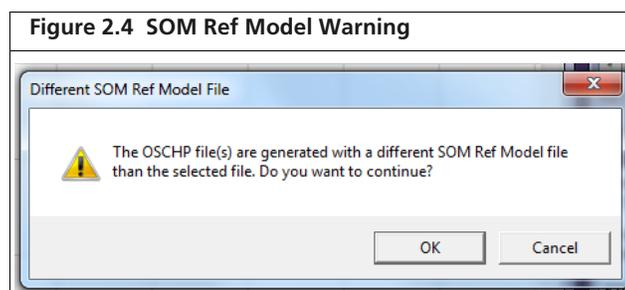
In most cases, the Viewer automatically chooses the appropriate location of the Reference Model file needed, as it is based on the OSCHP files you selected.

2. Click to select the Reference Model file, then click **Open**.

If your Reference Model File is not listed in the initial window, use the File window to navigate to the required Reference Model file, then click **Open**.

The Ref Model field is now populated.

A warning may appear if the selected SOM Ref Model is not the same as the one used to create the OSCHP files. (Figure 2.4)



- Click **Yes** to continue.
- Click **No** to remove the current SOM Ref Model File.

Assigning an Annotation File

! **IMPORTANT:** The annotation file provides information about each marker. Without this file, only the Probeset Name is available.

To assign the Annotation file:

1. Click **Select**.

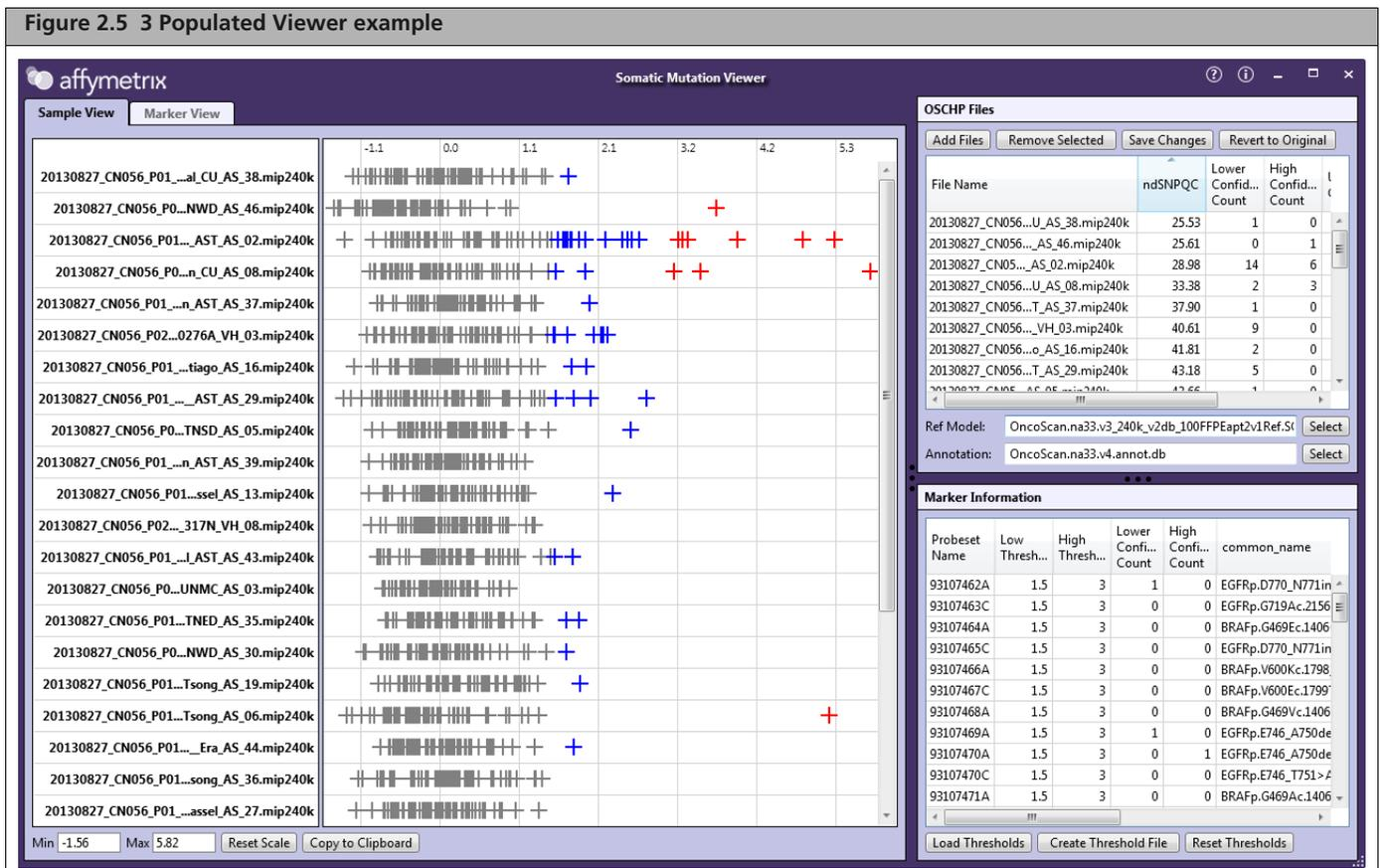
A File window appears.

In most cases, the Viewer automatically chooses the appropriate location of the Annotation file needed, as it is based on the OSCHP files you selected.

2. Click to select the Annotation file, then click **Open**.

If your Annotation File is not listed in the initial window, use the File window to navigate to the required Annotation file, then click **Open**.

The Somatic Mutation Viewer's 3 window panes are now fully populated. (Figure 2.5)



Using the Viewer

Different OSCHP files may have used different thresholds and/or somatic mutation reference model files. For display purposes, Somatic Mutation Viewer loads the thresholds from the first OSCHP file loaded.

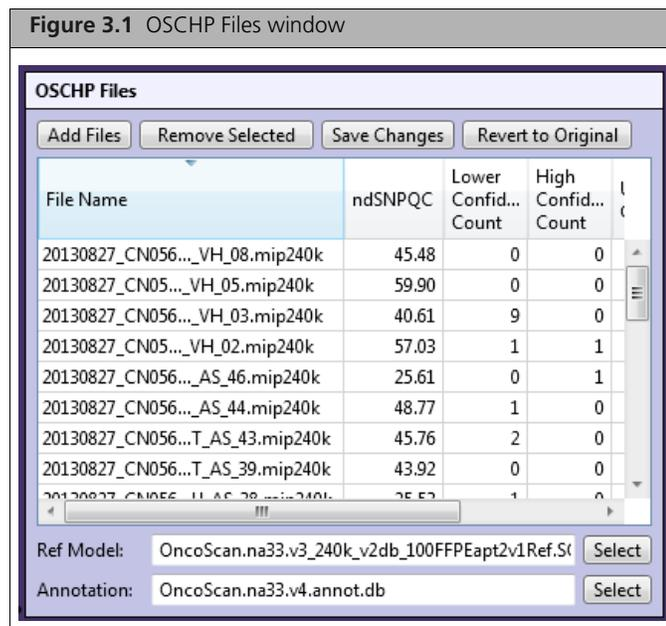
It is important that all OSCHPs must originate from the same SOM RefModel. If the thresholds are different, a message appears stating the first OSCHP thresholds will be used.

Poorer quality samples will have more false positive calls. Affymetrix recommends that you exclude samples that do not meet recommended QC thresholds, and any additional under-performing samples, from the visualizations and further analysis.

In the OSCHP Files table, sort on ndSNPQC to sort the samples according to their quality. Examine these samples in the Sample View. Consider removing the samples from the study that appear to have a substantial number of false positive calls.

OSCHP Files Window

The OSCHP Files window displays your OncoScan Console generated OSCHP files, the Reference Model File, and Annotation file. (Figure 3.1)



OSCHP File Window Columns

The Marker Information window (Figure 3.1) displays the following columns (from left to right):

- **File Name:** Displays the OSCHP file name.

- **ndSNPQC (SNP Quality Control of Normal Diploid Markers):** The metric SNPQC is a measure of how well genotype alleles are resolved in the microarray data. ndSNPQC is the same metric but only applied to normal diploid markers (that is those that have been determined to have Copy Number=2 in the sample). Larger ndSNPQC values are better.
- **Lower Confidence Count:** In the OSCHP Files table, this is the count of ProbeSets for the OSCHP that have a MutCall reporting “Lower confidence,” describing the likelihood that the mutation is present.
- **High Confidence Count:** In the OSCHP Files table, this is the count of ProbeSets for the OSCHP that have a MutCall reporting “High confidence,” describing the likelihood that the mutation is present.
- **Undetected Count:** This is the count of ProbeSets for the OSCHP that have a MutCall reporting “Undetected,” describing the likelihood that the mutation is not present.
- **Error Message:** Reports Somatic Mutation Viewer errors associated with the OSCHP file.

Adding and Removing OSCHP Files

To add OSCHP files to this window:

1. Click **Add Files**.
A File window appears.
2. Search within the File window that appears by default, or navigate to another folder location. Single click, Ctrl click, or Shift click (to select multiple OSCHP files).
3. Click **Open**.
The additional OSCHP files are now added.

To remove OSCHP files from this window:

1. Single click, Ctrl click, or Shift click (to select multiple OSCHP files).
2. Click **Remove Selected**.
The file(s) are removed.

Sorting OSCHP File Window Columns

To sort a column:

1. Click on a header.
The column is now sorted in an ascending order.
2. Click on the header again to reverse the sorting order.

To move a column:

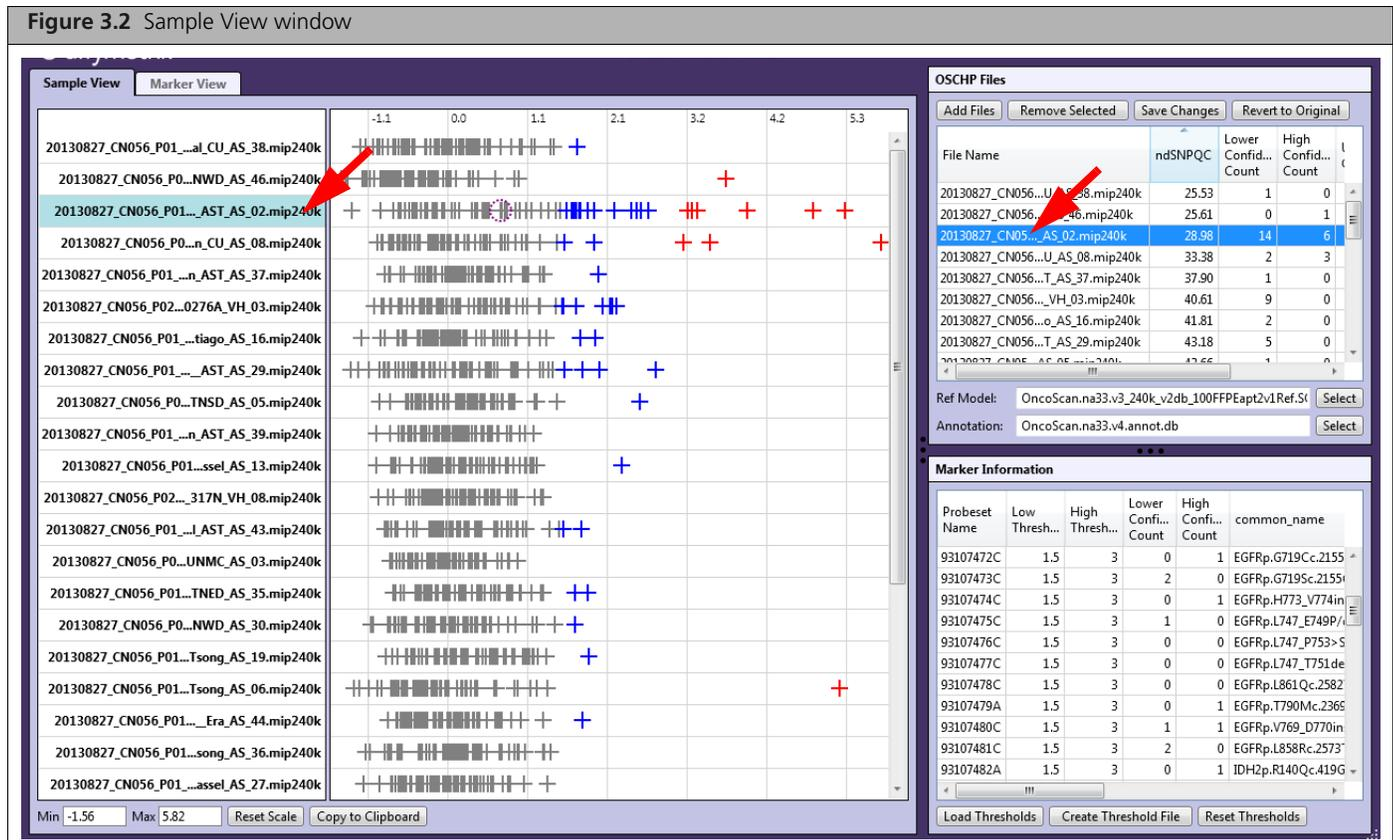
1. Click on a header, then drag it to a desired position within the OSCHP Files window.
2. Release the mouse button.
The header now resides at its new location within the OSCHP Files window.

Sample View Tab Window



TIP: The OSCHP Files window works in sync with the Sample View pane. Clicking on a sample file name also highlights that file within the Sample View pane, as shown in [Figure 3.2](#).

Your data is also in sync with the Marker Information window. If you click on a marker in the Sample View, that marker is highlighted in the Marker Information window, as shown in [Figure 3.2](#).



Using the Sample View Tab Window

The Sample View is **read-only**.

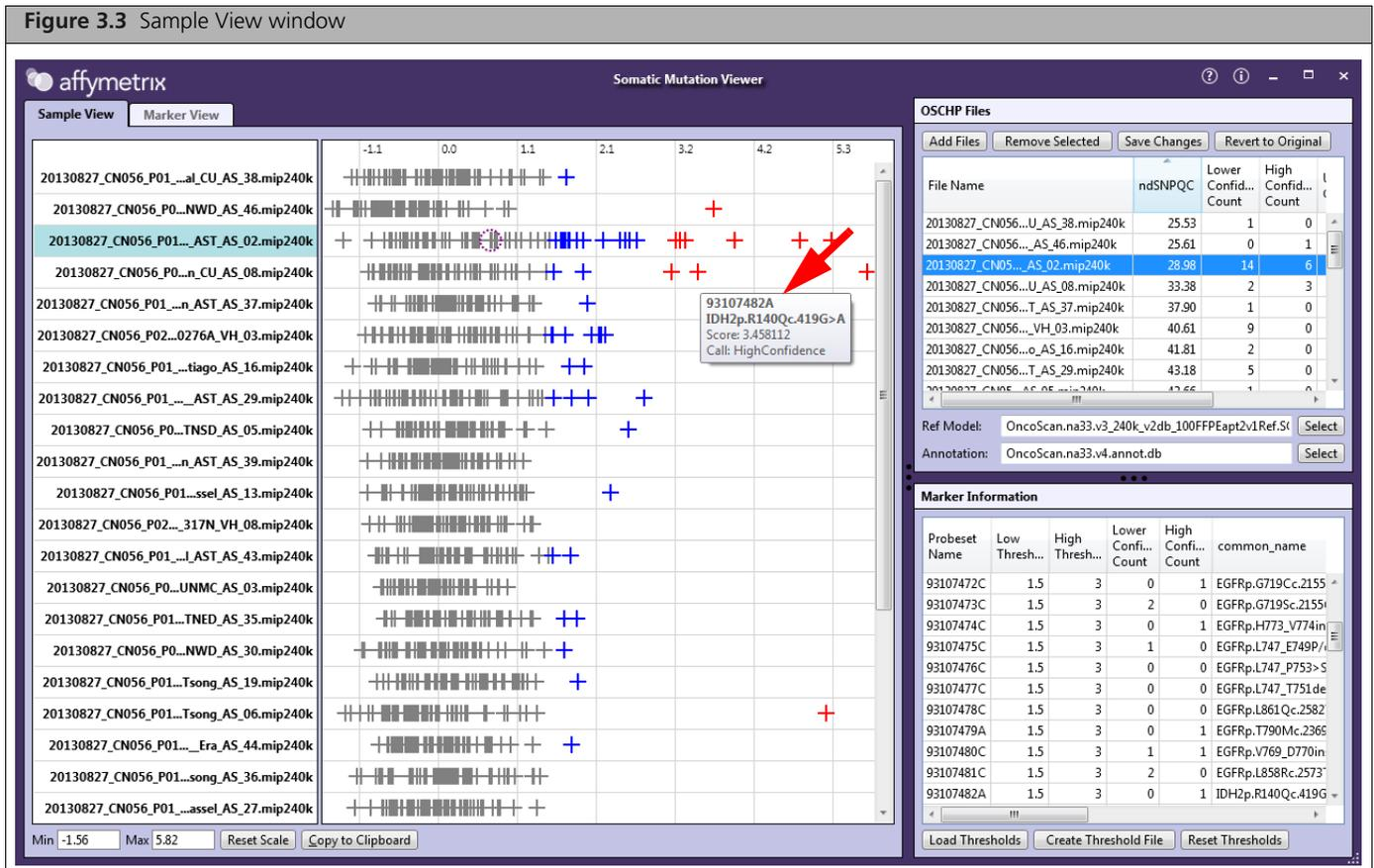
To view the Calls from an OSCHP File:

1. Click on the desired OSCHP file from the OSCHP File window pane or from the left pane of the Sample View window, as shown in [Figure 3.2](#).

To view specific calls:

1. Mouse over a call to reveal its score and Call property. Click on the call to see its Probeset Name and properties within the Marker Information window, as shown in [Figure 3.3](#).

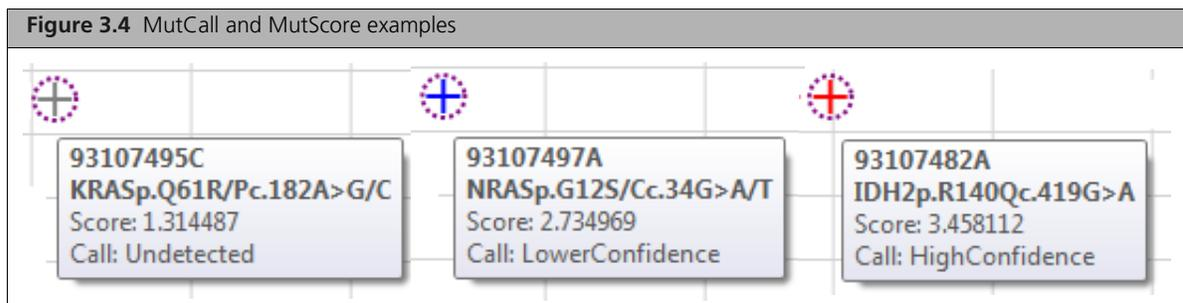
Figure 3.3 Sample View window



MutScores (Scores) and MutCalls (Calls) - Overview

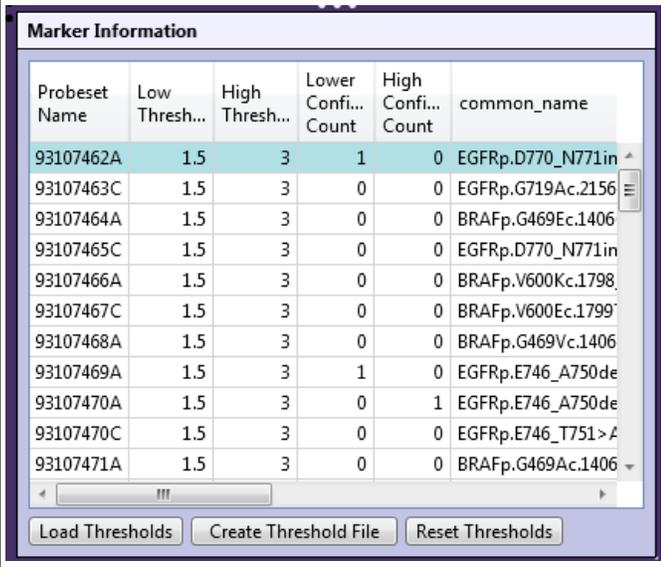
The x-axis signal is the MutScore from the OSCHP file. A MutScore is a measure of the signal response of the marker relative to the expected signal distribution of this marker in the absence of the mutation. It is calculated as follows:

- $\text{MutScore} = (\text{measured quantile normalized signal} - \text{median signal for this marker in the reference model file}) / (\text{95th percentile signal for this marker in the reference model file} - \text{median signal for this marker in the reference model file})$.
- A higher MutScore for the same marker indicates greater confidence that the mutation is present, and is correlated with higher % mutant allele. However, as each marker's signal is normalized to its own marker's reference signal distribution, it is not appropriate to compare MutScores between different markers to assign relative % mutant loads. Different markers will have different detection sensitivities.
- A MutCall is displayed as **Undetected** if the MutScore is below the Low Confidence threshold.
- A MutCall is reported as **HighConfidence** if greater than or equal to the High Confidence threshold.
- If the MutCall is equal to or greater than the Low Confidence threshold and is less than the High Confidence threshold, the MutCall is reported as **LowerConfidence**. (Figure 3.4)



Marker Information Window

Figure 3.5 Marker Information window.



Probeset Name	Low Thresh...	High Thresh...	Lower Confi... Count	High Confi... Count	common_name
93107462A	1.5	3	1	0	EGFRp.D770_N771in
93107463C	1.5	3	0	0	EGFRp.G719Ac.2156
93107464A	1.5	3	0	0	BRAFp.G469Ec.1406
93107465C	1.5	3	0	0	EGFRp.D770_N771in
93107466A	1.5	3	0	0	BRAFp.V600Kc.1798
93107467C	1.5	3	0	0	BRAFp.V600Ec.1799
93107468A	1.5	3	0	0	BRAFp.G469Vc.1406
93107469A	1.5	3	1	0	EGFRp.E746_A750de
93107470A	1.5	3	0	1	EGFRp.E746_A750de
93107470C	1.5	3	0	0	EGFRp.E746_T751>A
93107471A	1.5	3	0	0	BRAFp.G469Ac.1406

Buttons: Load Thresholds, Create Threshold File, Reset Thresholds

The Marker Information window (Figure 3.5) displays the following columns (from left to right):

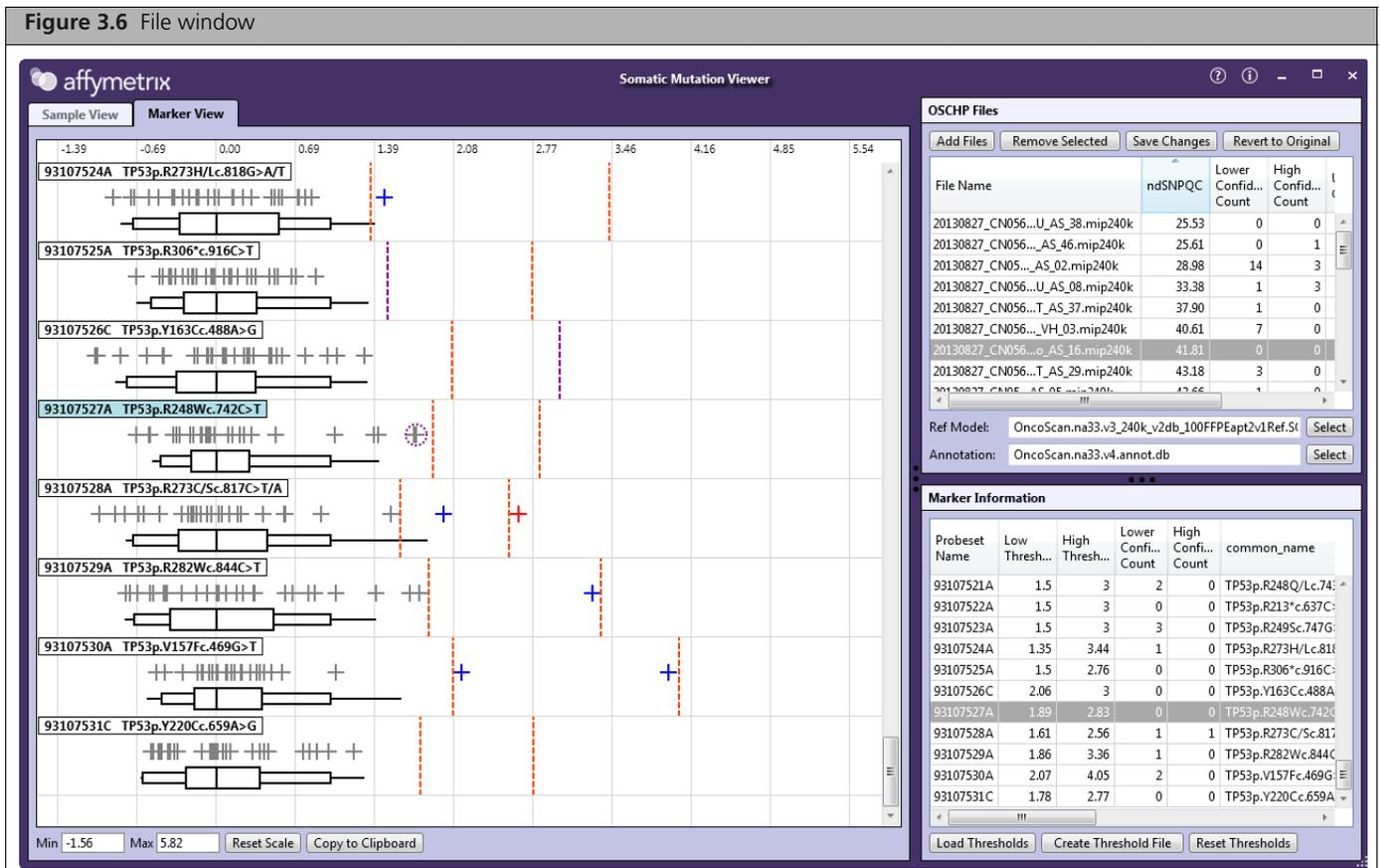
- **Probeset Name:** Affymetrix identifier for the marker.
- **Low Threshold:** Lower confidence MutScore threshold. Measurements with a MutScore below this value are called “Undetected”. Measurements equal to or greater than this threshold but less than the High Threshold are called “Lower confidence,” describing the likelihood that the mutation is present.
- **High Threshold:** High confidence MutScore threshold. Measurements equal to or greater than this threshold are called “High confidence,” describing the likelihood that the mutation is present.
- **Lower Confidence Count:** In the Marker Information table, this is the count of OSCHP files for the ProbeSet that have a MutCall reporting “Lower confidence.”
- **High Confidence Count:** In the Marker Information table, this is the count of OSCHP files for the ProbeSet that have a MutCall reporting “High confidence.”
- **common_name:** Abbreviated description of the mutations to which this ProbeSet is known to respond. The name has the form [Gene]:[amino acid change for mutation]:[cDNA change for mutation]. In the event that the ProbeSet cannot differentiate among multiple mutations to which it can respond, the slash (/) delimits the multiple known mutations.
- **chr_ID:** Chromosome of the mutation.
- **start:** Starting genomic position of the mutation.
- **stop:** Ending genomic position of the mutation.

- **cosmic_id:** The identifier of the mutation as listed in the COSMIC database, which is a catalogue of somatic mutations in cancer. More information on these mutations can be found at <http://cancer.sanger.ac.uk>
- **channel:** The CEL file from which the signal is measured. "A" is the AT CEL, "C" is the GC CEL.



TIP: The Marker Information window works in sync with the Marker View pane. Clicking on a sample file name also highlights that file within the Marker View pane, as shown in [Figure 3.6](#).

Figure 3.6 File window



Editing Thresholds in the Marker Information Window

To edit the Low and/or High Thresholds:

1. Click to highlight the Probeset marker you want to edit.

- Click on the current *Low Threshold* value field.

Figure 3.7 Changing Low Threshold value example 1

Marker Information						
Probeset Name	Low Thresh...	High Thresh...	Lower Confid... Count	High Confid... Count	chr_id	start
93107469A	1.5	3	1	0	7	55242464
93107470A	1.5	3	0	1	7	55242465
93107470C	1.5	3	0	0	7	55242466

- Enter a new *Low Threshold* value.

Figure 3.8 Changing Low Threshold value example 2

Marker Information						
Probeset Name	Low Thresh...	High Thresh...	Lower Confid... Count	High Confid... Count	chr_id	start
93107469A	1.5	3	1	0	7	55242464
93107470A	1.8	3	0	1	7	55242465
93107470C	1.5	3	0	0	7	55242466

- Press **Enter**.
The Low Threshold field value is now changed. This threshold value change (represented by vertical dotted lines) is also reflected in the Marker View window.
- If you want to change the current *High Threshold* value field, click on the current *High Threshold* value field.
- Enter a new *High Threshold* value.
- Press **Enter**.
The High Threshold field value is now changed. This threshold value change (represented by vertical dotted lines) is also reflected in the Marker View window.

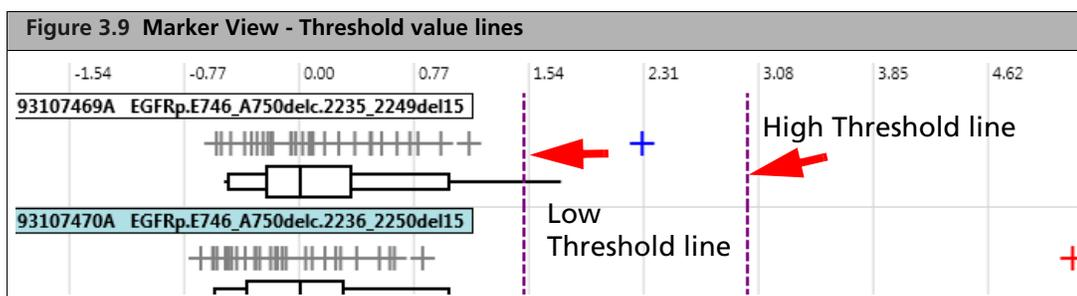
Marker View Tab Window

Thresholds for individual markers can be changed either by dragging the vertical dashed threshold lines in the Marker View, or by manually editing the Low and High Threshold values in the Marker Info table.

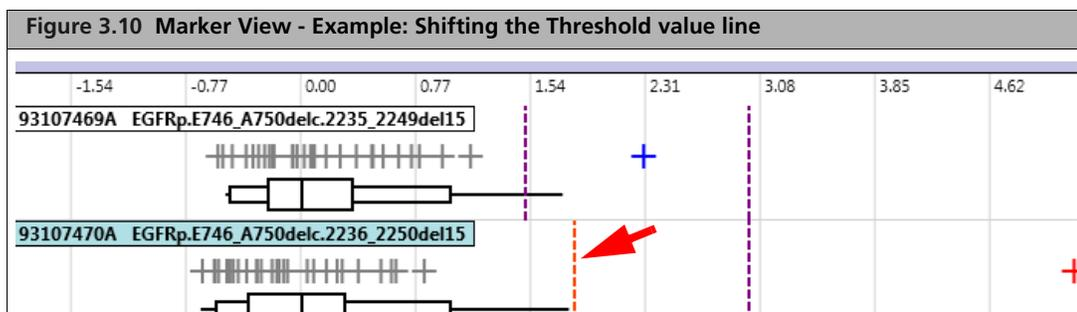
Editing Thresholds in the Marker View Tab Window

To edit the Low and/or High Thresholds:

1. Click to highlight the Probeset marker you want to edit.
Low and High Thresholds are represented by vertical dotted lines, as shown in [Figure 3.9](#).



2. Click on the current *Low Threshold* line, drag the line to your desired location, then release the mouse button. ([Figure 3.10](#)).



The Low Threshold field value is now changed. This threshold value change is also reflected in the Marker Information window's corresponding Low Threshold value field.

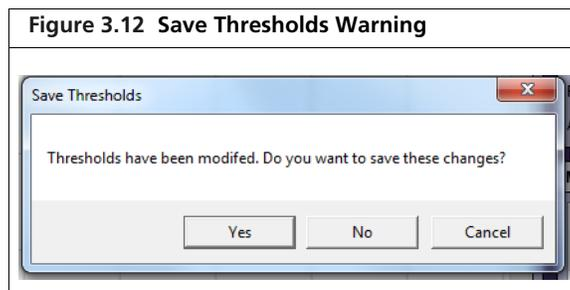
3. If you want to change the current *High Threshold* value field, click on the current *High Threshold* line, drag the line to your desired location, then release the mouse button.
The High Threshold field value is now changed. This threshold change is also reflected in the Marker Information window's corresponding High Threshold value field.

A warning may appear when adding or removing an OSCHP file (after its thresholds have been changed). (Figure 3.11)



Do one of the following:

- Click **OK** to continue.
- Clicking **Cancel** prompts the following message: (Figure 3.12)



Do one of the following:

1. Click: **Yes** - Saves the updated calls and thresholds to the OSCHP file and closes the viewer.
2. Click **No** - Does not save the updated thresholds and any changes to the calls and closes the viewer.
3. Click **Cancel** - Does not close the Viewer.

Saving your New Thresholds



TIP: The Viewer always retains your original thresholds. At any time, click *Reset Thresholds* to restore your original threshold values.

To save your new thresholds:

1. At the Marker Information window, click **Create Thresholds**.
A File window appears.
2. Navigate to a desired save location, enter a filename, then click **Save**.
Your new Thresholds file is saved as a Tab-delimited text file. You can then load this Thresholds file into Somatic Mutation Viewer at another time, or you can load it into OncoScan Console when setting up additional analyses of CEL files.

To load thresholds:

1. At the Marker Information window, click **Load Thresholds**.

A File window appears.

2. Navigate to the Threshold's location, click on its filename, then click **Open**. Your Thresholds file's properties appear in the Marker Information window.

Viewing Tools

- For optimum viewing, each Somatic Mutation Viewer window pane can be easily resized
- The Marker and Sample Views feature a taskbar that can customize your desired view even further.
- The Marker View features a Box Whisker Plot graphic for each Probeset.

Resizing Window Panes

To resize a window pane:

1. Click on the edge of a window you want to resize, then drag it to the size you want.

Taskbar Options

The Marker and Sample Views feature a taskbar (bottom left) that enables you to do the following: (Figure 3.13)



- **Min:** Enter your minimum (starting point) scale size of the currently displayed data.
- **Max:** Enter your maximum (finishing point) scale size of the currently displayed data.
- **Reset Scale:** Returns your view to its original scaled state - as it was - the first time the files were loaded into the Viewer.
- **Copy to Clipboard:** Click to save the current (Marker or Sample) view to your Clipboard for pasting in another application (as a .PNG file).

Box Whisker Plot Graphic

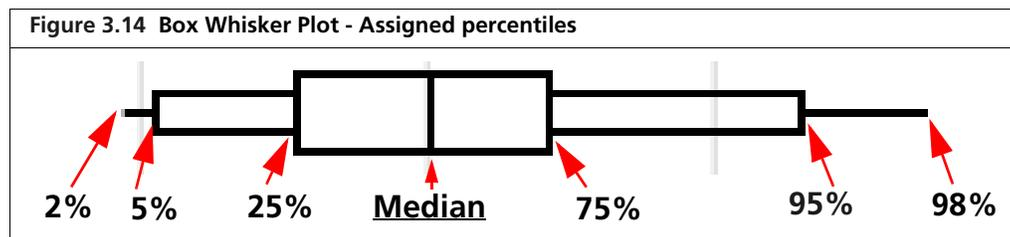
If you have selected a reference model file, the reference signal distribution for each marker in the absence of any mutation is displayed as a box plot in the Marker View.

The percentile values marked by the box plot represent the 2%, 5%, 25%, 50%, 75%, 95%, and 98% percentiles of the signal distribution of this marker in the reference. The 50%ile value of the reference will have a MutScore = 0, and the 95%ile value will have a MutScore = 1.

Make sure the reference model file you select is the one used to generate the OSCHP file results you are displaying.

Consider utilizing this plot graphic when determining placement of new threshold values.

Even though each Probeset's plot is uniquely sized, its assigned percentile locations are constant, as shown in [Figure 3.14](#).



Saving your Edited OSCHP Files



TIP: When you use Somatic Mutation Viewer to save changed thresholds and calls to your OSCHP files, the original values are retained in the OSCHP files.

At any time, click *Revert to Original* to restore your original thresholds and calls.

To save your edited OSCHP file set:

1. At the OSCHP Files window, click **Save Changes**.
A green progress bar appears and your newly edited OSCHP file set is now saved.

To restore your original OSCHP file set:

1. At the OSCHP Files window, click **Revert to Original**.
A green progress bar appears and your original OSCHP file set is now restored.