

A large, metallic, silver-colored DNA double helix structure that spirals diagonally from the bottom left towards the top right of the page.

## *User Manual*

---

# Axiom CNV Summary Tool

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

## Trademarks

Affymetrix®, Axiom®, GeneChip®, NetAffx®, Command Console®, Powered by Affymetrix™, GeneChip-compatible™, Genotyping Console™, DMET™, GeneTitan®, 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.

---

<b>Chapter 1</b>	<b>Introduction . . . . .</b>	<b>2</b>
	Prerequisites . . . . .	2
	Recommended System Requirements . . . . .	2
	Installing Axiom CNV Summary Tool . . . . .	2
	Starting Axiom CNV Summary Tool . . . . .	2
	Overview of Workflow and Useful Tips . . . . .	4
	Overview of Workflow . . . . .	4
	Useful Tips . . . . .	5

---

<b>Chapter 2</b>	<b>Axiom CNV Summary Tool . . . . .</b>	<b>7</b>
	GTC Users . . . . .	7
	APT Users . . . . .	7
	Using the Axiom CNV Summary Tool . . . . .	8
	Select Gender File. . . . .	9
	Reference File (Optional) . . . . .	9
	Running the Axiom CNV Summary Tool . . . . .	10
	Retrieving the Axiom CNV Summary Tool Data . . . . .	10
	Ways to Use the Axiom CNV Summary Tool Data . . . . .	10
	Subsequent Analyses . . . . .	10
	Viewing Data in the Axiom CNV Viewer . . . . .	11
	Further Copy Number Analysis Using BioDiscovery's Nexus Software . . . . .	16

---

<b>Chapter 3</b>	<b>Performing GC Correction in Nexus . . . . .</b>	<b>18</b>
------------------	----------------------------------------------------	-----------

---

<b>Appendix A</b>	<b>Calculations of log<sub>2</sub> Ratios and B allele Frequencies . . . . .</b>	<b>23</b>
	Computing log <sub>2</sub> Ratio . . . . .	23
	B allele Frequencies (BAF) . . . . .	24

## Introduction

This document provides the general guidelines on how to use the **Axiom CNV Summary Tool**.

The **Axiom CNV Summary Tool** generates input files for BioDiscovery Nexus using Axiom data.

The included **Axiom CNV Viewer** allows you to view the data generated by the Axiom CNV Summary Tool. To use the Viewer, see [Viewing Data in the Axiom CNV Viewer on page 11](#)

## Prerequisites

Before you can use the Axiom CNV Tools you must have:

- An appropriate Annotation File (\*annot.db) downloaded from the Affymetrix NetAffx website. (User Account and Password required)
- Good quality genotyping data processed through the Affymetrix Genotyping Console™ (GTC) or Affymetrix Power Tools™ (APT) application.

## Recommended System Requirements

64-bit Operating System	Speed	Memory (RAM)	Available Disk Space	Web Browser
Microsoft Windows 7 Professional	2.83 GHz Pentium Quad Core Processor	16 GB	150 GB HD + Data Storage	IE 8.0 and above
Microsoft Windows 8 Professional	2.83 GHz Pentium Quad Core Processor	16 GB	150 GB HD + Data Storage	IE 8.0 and above

## Installing Axiom CNV Summary Tool

To install the Axiom CNV Summary Tool:

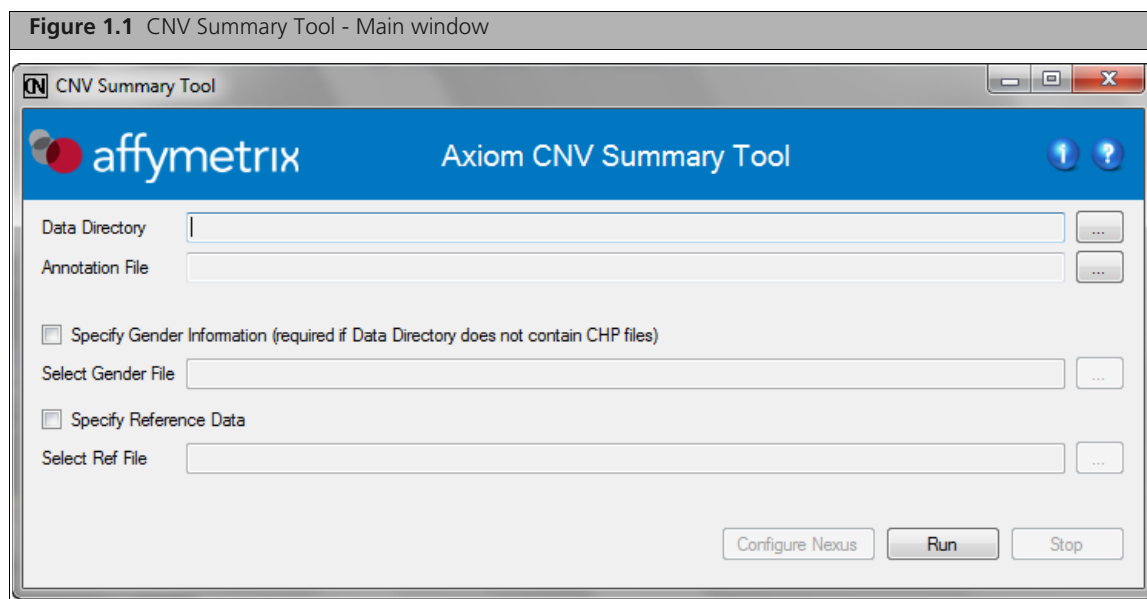
1. Go to **www.affymetrix.com** and navigate to the following location:  
**Home > Products > Microarray Solutions > Instruments and Software > Software >**
2. Locate and download the zipped **Axiom CNV Summary Tool** software package.
3. Unzip the file, then double-click **AxiomCNVSetup.exe** to install it.
4. Follow the installer's instructions.

## Starting Axiom CNV Summary Tool

To start the CNV Summary Tool:

1. Click **Start -> All Programs -> Affymetrix -> Axiom CNV Tools**.

2. Locate and click on **Axiom CNV Summary Tool**.  
The application opens. (Figure 1.1)

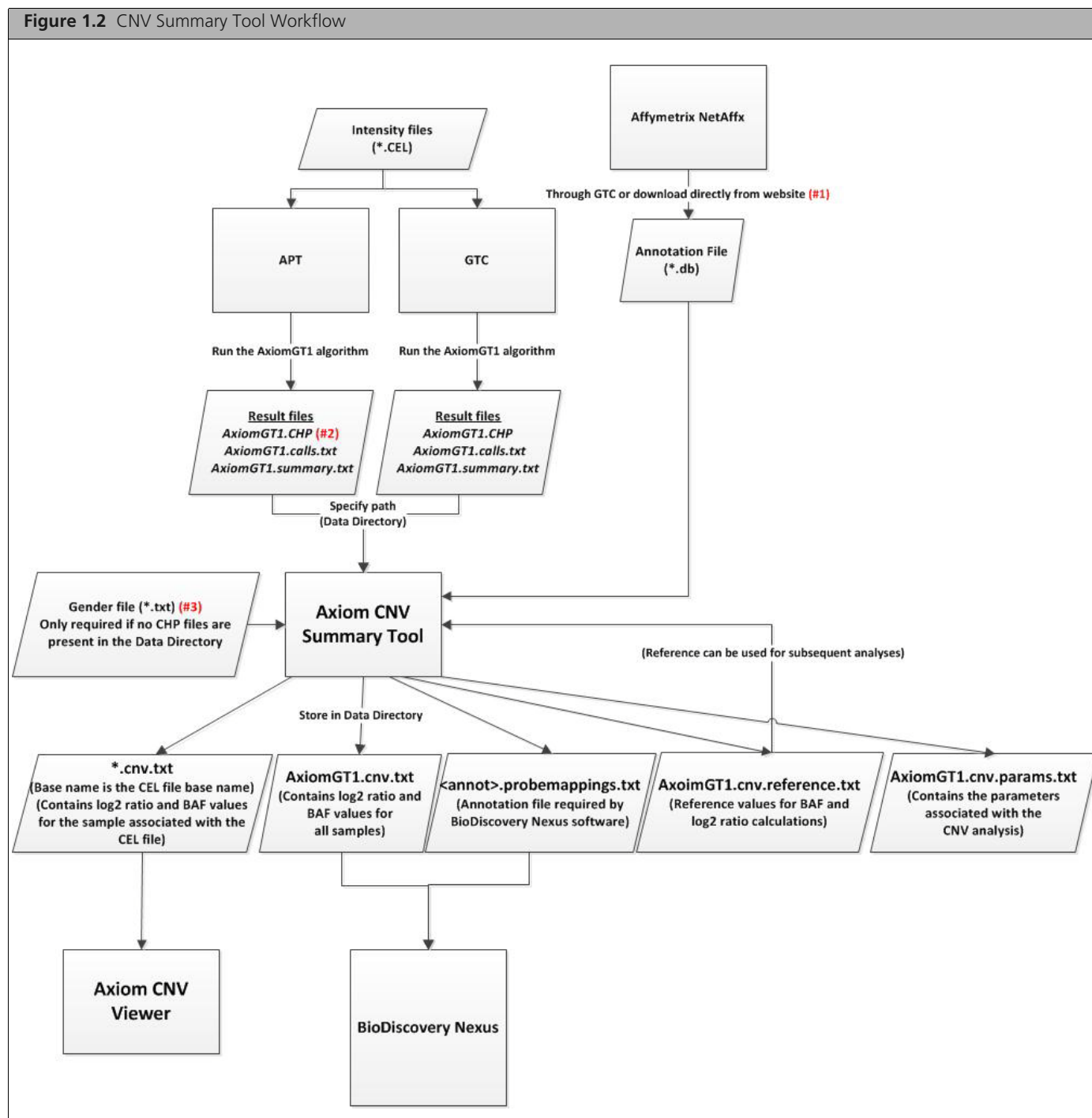


## Overview of Workflow and Useful Tips

### Overview of Workflow

Please review the diagram below. It shows the data analysis workflow starting with GTC or APT and ending with data for use with BioDiscovery Nexus.

The file types that are required and generated by the Axiom CNV Summary Tool are also noted in this diagram. (Figure 1.2)



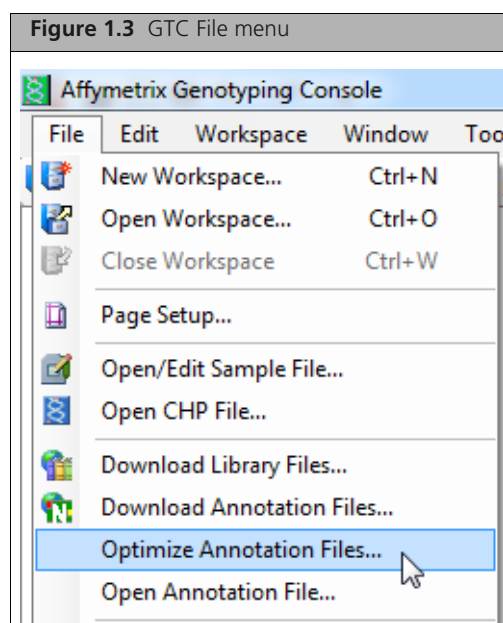
## Useful Tips

### Tip #1

When you download an annotation file using GTC, the annotation file is automatically indexed for optimum processing.

If your annotation file was not downloaded from GTC, you can optimize it for faster processing using the GTC software, as follows:

Click **File** -> **Optimize Annotation File...** (Figure 1.3)



### Tip #2

To create CHP files in ATP, use the parameter: **--cc-chp-output**

### Tip #3

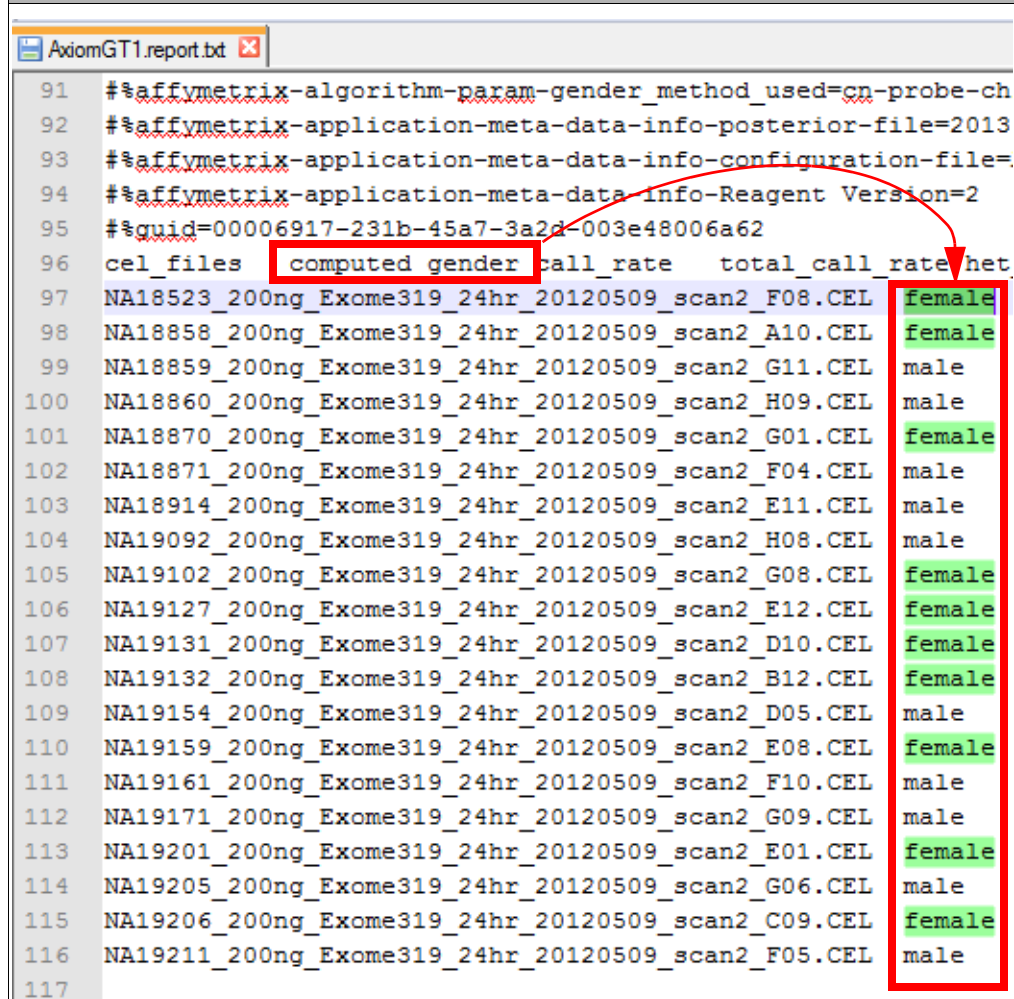
You can create your own gender.txt file by using the information in the *AxiomGT1.report.txt* file. This file is located in your *Genotyping Results* folder.

The second column in the report.txt file is *computed\_gender*. (Figure 1.4)

Select, then copy the gender information you want to use from the report.txt file, then paste it in a new *gender.txt* file.

See [Figure 2 on page 9](#) for gender.txt file format requirements, then follow steps; [To include gender values for all samples: on page 9](#).

Figure 1.4 Axiom GT1.report.txt file example



```

91  #%affymetrix-algorithm-param-gender_method_used=cn-probe-ch
92  #%affymetrix-application-meta-data-info-posterior-file=2013
93  #%affymetrix-application-meta-data-info-configuration-file=
94  #%affymetrix-application-meta-data-info-Reagent Version=2
95  #%guid=00006917-231b-45a7-3a2d-003e48006a62
96  cel_files  computed gender  call_rate  total_call_rate  net
97  NA18523_200ng_Exome319_24hr_20120509_scan2_F08.CEL  female
98  NA18858_200ng_Exome319_24hr_20120509_scan2_A10.CEL  female
99  NA18859_200ng_Exome319_24hr_20120509_scan2_G11.CEL  male
100 NA18860_200ng_Exome319_24hr_20120509_scan2_H09.CEL  male
101 NA18870_200ng_Exome319_24hr_20120509_scan2_G01.CEL  female
102 NA18871_200ng_Exome319_24hr_20120509_scan2_F04.CEL  male
103 NA18914_200ng_Exome319_24hr_20120509_scan2_E11.CEL  male
104 NA19092_200ng_Exome319_24hr_20120509_scan2_H08.CEL  male
105 NA19102_200ng_Exome319_24hr_20120509_scan2_G08.CEL  female
106 NA19127_200ng_Exome319_24hr_20120509_scan2_E12.CEL  female
107 NA19131_200ng_Exome319_24hr_20120509_scan2_D10.CEL  female
108 NA19132_200ng_Exome319_24hr_20120509_scan2_B12.CEL  female
109 NA19154_200ng_Exome319_24hr_20120509_scan2_D05.CEL  male
110 NA19159_200ng_Exome319_24hr_20120509_scan2_E08.CEL  female
111 NA19161_200ng_Exome319_24hr_20120509_scan2_F10.CEL  male
112 NA19171_200ng_Exome319_24hr_20120509_scan2_G09.CEL  male
113 NA19201_200ng_Exome319_24hr_20120509_scan2_E01.CEL  female
114 NA19205_200ng_Exome319_24hr_20120509_scan2_G06.CEL  male
115 NA19206_200ng_Exome319_24hr_20120509_scan2_C09.CEL  female
116 NA19211_200ng_Exome319_24hr_20120509_scan2_F05.CEL  male
117

```



# Axiom CNV Summary Tool

Before you can run the Axiom CNV Summary Tool, you must first process the CEL files through one of the following software applications:

- Affymetrix Genotyping Console (GTC) v4.1.4 (or higher)
- Affymetrix Power Tools (APT)

## GTC Users

Do the following:

1. Run the AxiomGT1 algorithm through GTC (using your \*.CEL Intensity files) as you normally would.

GTC produces the following Result files:

- CHP files
- AxiomGT1.calls.txt
- AxiomGT1.summary.txt

## APT Users

Do the following:

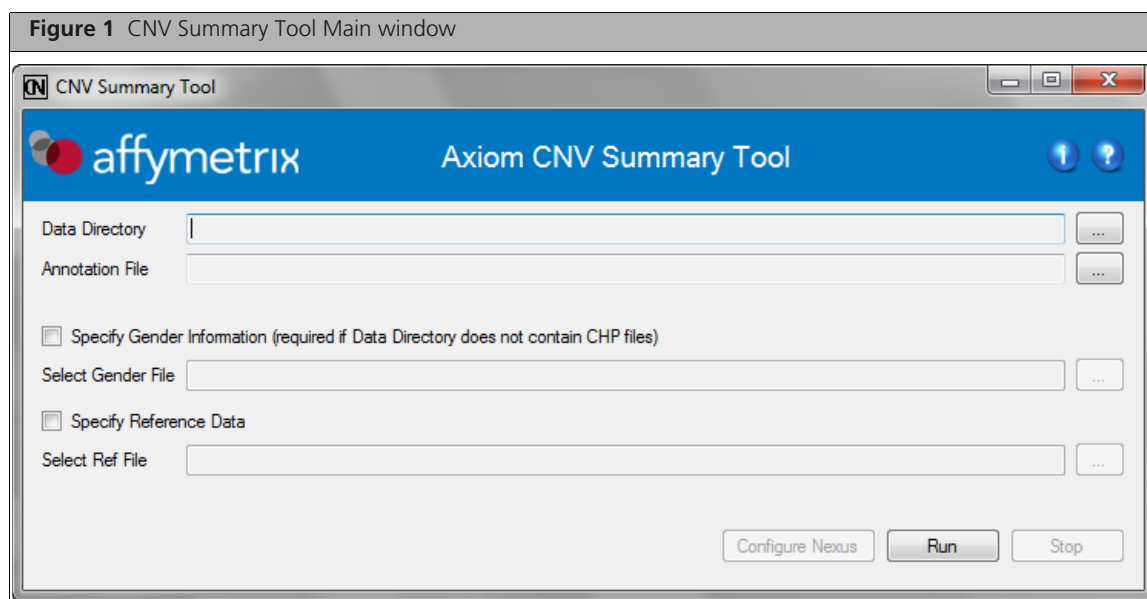
1. Run the AxiomGT1 algorithm through APT (using your \*.CEL Intensity files) as you normally would.

APT produces the following Result files:

- AxiomGT1.calls.txt
- AxiomGT1.summary.txt
- CHP files (APT's **--cc-chp-output** option must be enabled)

## Using the Axiom CNV Summary Tool

From the Axiom CNV Summary Tool window: (Figure 1)



1. Click the **Data Directory** Browse button.
2. Navigate to the folder that contains your input data files (\*AxiomGT1.CHP, AxiomGT1.calls.txt, and AxiomGT1.summary.txt files), then single click, Ctrl click, Shift click, or Ctrl-A (to select multiple files).
3. Click **OK**.  
The Data Directory path is now populated.
4. Click the **Annotation File** Browse button.
5. Navigate to the folder that contains the annotation file you downloaded earlier from [www.affymetrix.com](http://www.affymetrix.com).



**NOTE:** Annotation files are array specific. If you are running an analysis for a specific array, make sure you use the appropriate annotation file.

Annotation files for *Axiom myDesign* arrays are provided directly to you from Affymetrix (for each custom array designed).

When you download an annotation file using GTC, the annotation file is automatically indexed for optimum processing. If your annotation file was not downloaded from GTC, see [Tip #1 on page 5](#).

6. Click to select the annotation file, then click **OK**.  
The Annotation File path is now populated

## Select Gender File.

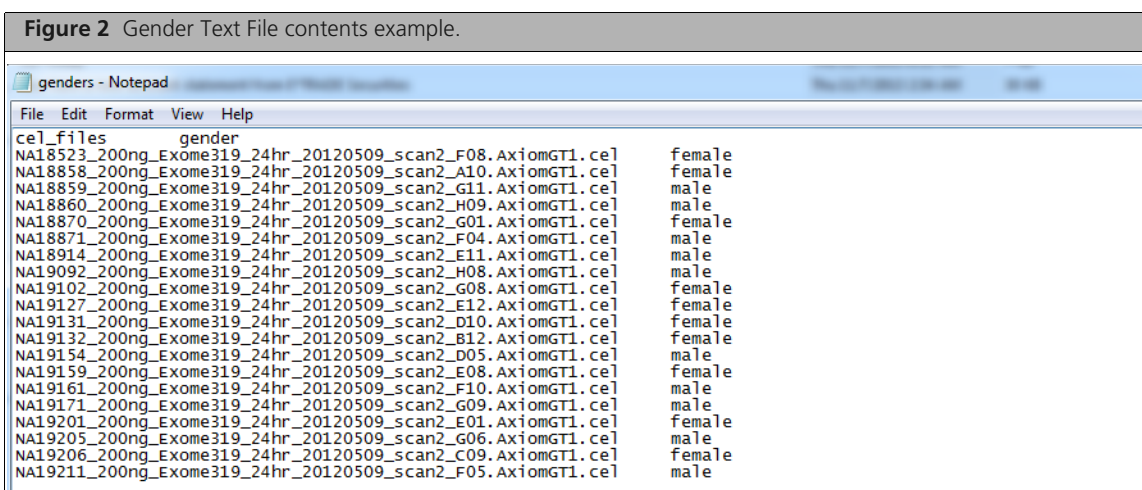
**▲ IMPORTANT:** If there are no CHP files present in your Data Directory folder, you **MUST** include gender values for all samples. If you are using CHP files, skip this *Select Gender File* option.

### To include gender values for all samples:

1. Click the **Specify Gender information** checkbox.
2. Click the **Gender File** Browse button.
3. Navigate to the folder that contains your Gender files.

**📄 NOTE:** Your gender file must be a tab-delimited text file with 2 columns. Its first column header must be *cel\_files*. The second column header must be *gender*, as shown in [Figure 2](#).

**Figure 2** Gender Text File contents example.



cel_files	gender
NA18523_200ng_Exome319_24hr_20120509_scan2_F08.AxiomGT1.cel	female
NA18858_200ng_Exome319_24hr_20120509_scan2_A10.AxiomGT1.cel	female
NA18859_200ng_Exome319_24hr_20120509_scan2_G11.AxiomGT1.cel	male
NA18860_200ng_Exome319_24hr_20120509_scan2_H09.AxiomGT1.cel	male
NA18870_200ng_Exome319_24hr_20120509_scan2_G01.AxiomGT1.cel	female
NA18871_200ng_Exome319_24hr_20120509_scan2_F04.AxiomGT1.cel	male
NA18914_200ng_Exome319_24hr_20120509_scan2_E11.AxiomGT1.cel	male
NA19092_200ng_Exome319_24hr_20120509_scan2_H08.AxiomGT1.cel	male
NA19102_200ng_Exome319_24hr_20120509_scan2_G08.AxiomGT1.cel	female
NA19127_200ng_Exome319_24hr_20120509_scan2_E12.AxiomGT1.cel	female
NA19131_200ng_Exome319_24hr_20120509_scan2_D10.AxiomGT1.cel	female
NA19132_200ng_Exome319_24hr_20120509_scan2_B12.AxiomGT1.cel	female
NA19154_200ng_Exome319_24hr_20120509_scan2_D05.AxiomGT1.cel	male
NA19159_200ng_Exome319_24hr_20120509_scan2_E08.AxiomGT1.cel	female
NA19161_200ng_Exome319_24hr_20120509_scan2_F10.AxiomGT1.cel	male
NA19171_200ng_Exome319_24hr_20120509_scan2_G09.AxiomGT1.cel	male
NA19201_200ng_Exome319_24hr_20120509_scan2_E01.AxiomGT1.cel	female
NA19205_200ng_Exome319_24hr_20120509_scan2_G06.AxiomGT1.cel	male
NA19206_200ng_Exome319_24hr_20120509_scan2_C09.AxiomGT1.cel	female
NA19211_200ng_Exome319_24hr_20120509_scan2_F05.AxiomGT1.cel	male

The Gender column (far right) ([Figure 2](#)) is not case-sensitive.

- For *Female* type: **Female, female, F, or 2**
- For *Male* type: **Male, male, M, or 1**
- To specify an *unknown* gender type: **unknown or 0**

If no gender was specified or the gender was specified other than the required naming conventions stated above, the gender entry will be treated as *unknown*.

4. Click to select the *gender.txt* file you want to use, then click **OK**.

The Gender File path is now populated.

## Reference File (Optional)

The choice of samples to be used as a reference is critical for accurate CNV detection because the log2ratio at a marker is computed by dividing the intensity of the marker by the median intensity of that marker in the chosen reference set, in log space. The reference set, therefore, should represent the normal copy number state for each marker. One approach is to create the reference based on the individuals genotyped on the plate, provided that for each marker the vast majority of individuals on the plate are expected to have normal copy number states. Another approach is to create a separate reference based

on individuals expected to have normal copy number states genotyped on different plates. If the latter approach is chosen the number of samples used for the reference should be as large as possible, preferably at least 100. The analysis can be carried out with any number of samples but will be less accurate for smaller reference sets.

#### Do the following to specify reference data:

1. Click the **Specify Reference Data** checkbox.
2. Click the **Select Ref File** Browse button.
3. Navigate to the folder that contains your reference data file.
4. Click to select the file you want, then click **OK**.

The Select Ref File path is now populated.

### Running the Axiom CNV Summary Tool

1. After your Axiom CNV Summary Tool data paths are set, click **Run**.

A green progress bar appears. Processing time varies depending on the amount of data you are processing, the number of SNPs on your array, and your system's hardware specifications.

After the data has been successfully processed, a message appears.

2. Click **OK** to acknowledge the message.

### Retrieving the Axiom CNV Summary Tool Data

The following files are produced and are stored in the **Data Directory** folder you assigned earlier:

- **\*.cnv.txt** - Base name is the CEL file base name. (This file contains log2 ratio and BAF values for the sample associated with the CEL file.)
- **AxiomGT1.cnv.txt** - Contains log2 ratio and BAF values for all samples.
- **<annot>.probemappings.txt** - Required by BioDiscovery's Nexus software.
- **AxiomGT1.cnv.reference.txt** - Reference values for BAF and log2 ratio calculations.
- **AxiomGT1.cnv.params.txt** - Contains the parameters associated with the CNV analysis.

## Ways to Use the Axiom CNV Summary Tool Data

### Subsequent Analyses

Use the newly generated **AxiomGT1.cnv.reference.txt** for additional analysis.

1. Click the **Specify Reference Data** checkbox.
2. Click the Select Ref File Browse button, navigate to your Data Directory folder, then click to select the file: **AxiomGT1.cnv.reference.txt**
3. After your Axiom CNV Summary Tool data paths are set, click **Run**.

A green progress bar appears. Allow time for your data to process.

After the data has been successfully processed, a message appears.

4. Click **OK** to acknowledge the message.

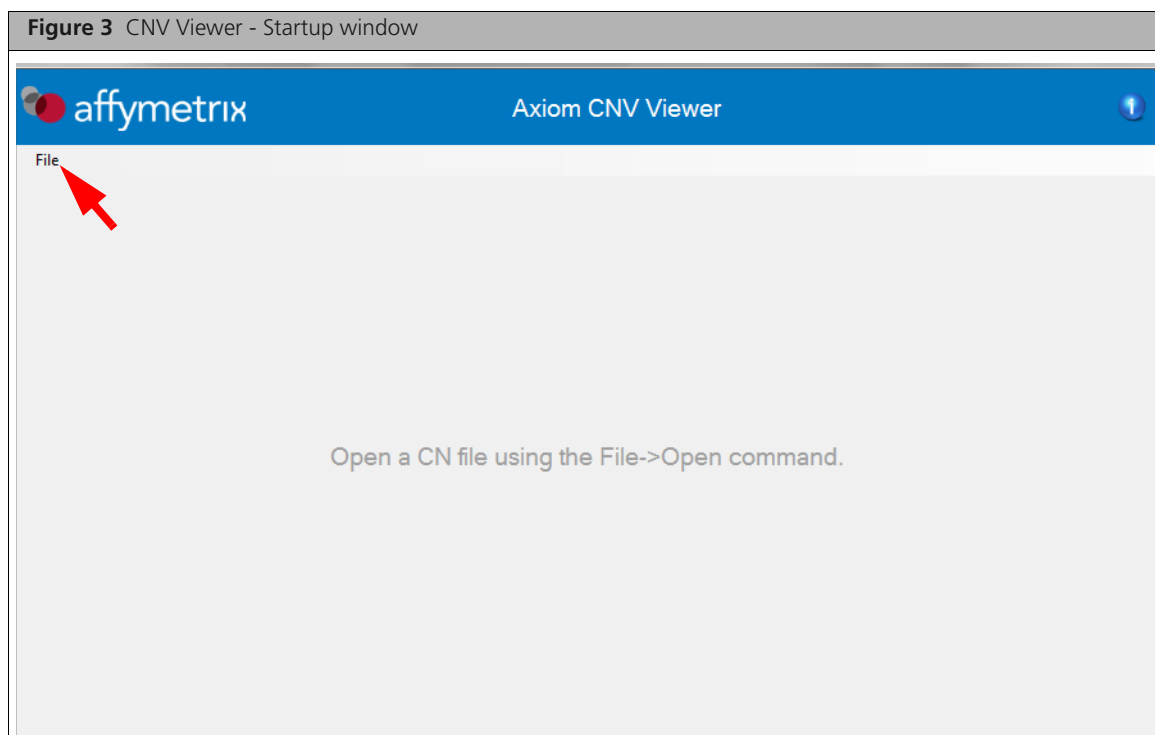
## Viewing Data in the Axiom CNV Viewer

View the newly generated **\*.cnv.txt** for additional analysis in the included Axiom CNV Viewer.

### To start the Axiom CNV Viewer:

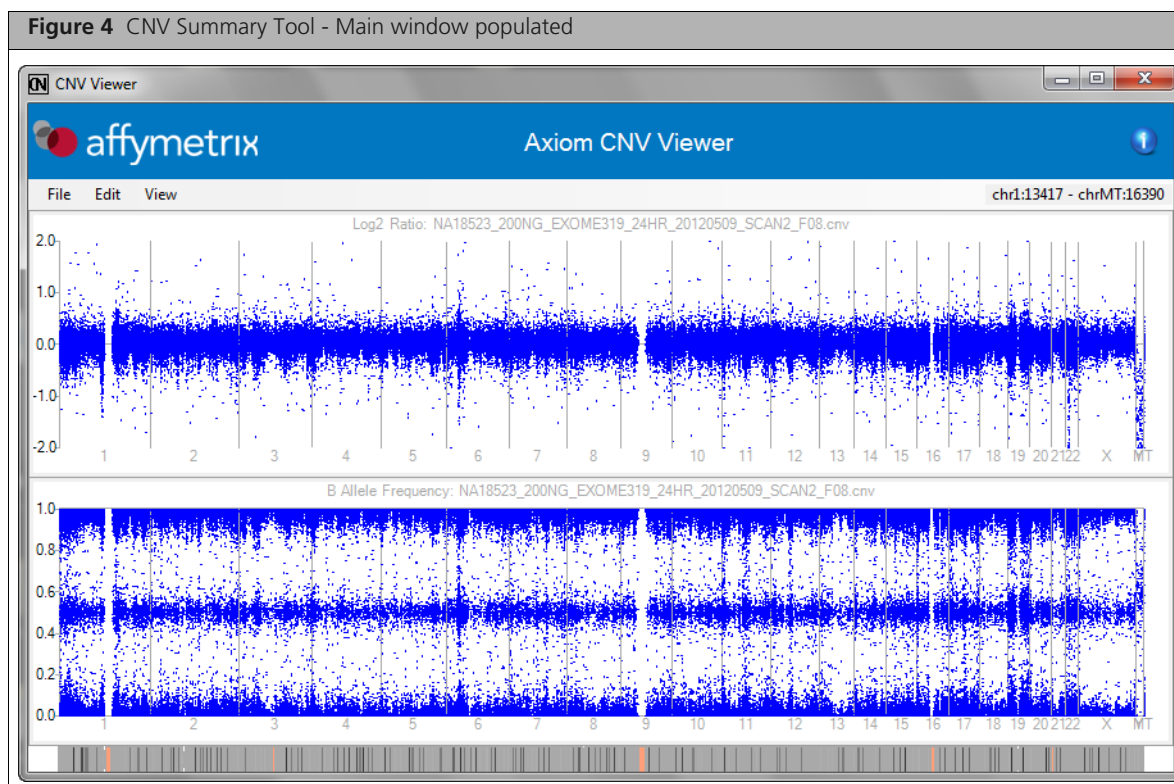
1. Click **Start** -> **All Programs** -> **Affymetrix** -> **Axiom CNV Tools**.
2. Locate and click on **Axiom CNV Viewer**.

The application opens. (Figure 3)



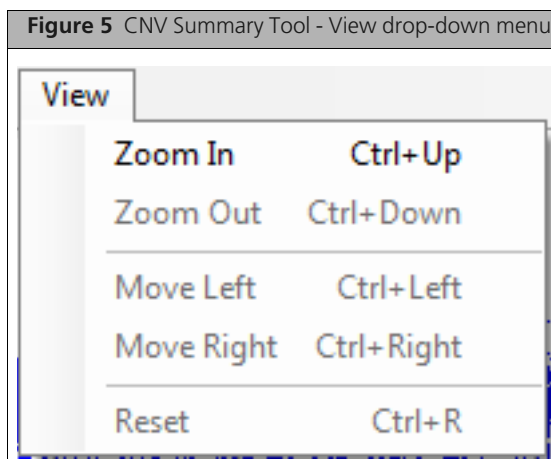
3. Click **File** -> **Open**.
4. Navigate to your Data Directory folder, then select the **\*.cnv.txt** file(s) you want to view.
5. Click **OK**.

The Viewer displays your data. (Figure 4)



**To customize the display view:**

1. Click **View**, then click to select one of the following viewing options or use the equivalent keyboard commands shown. (Figure 5)
2. Repeat the viewing command as needed to reach the desired view.



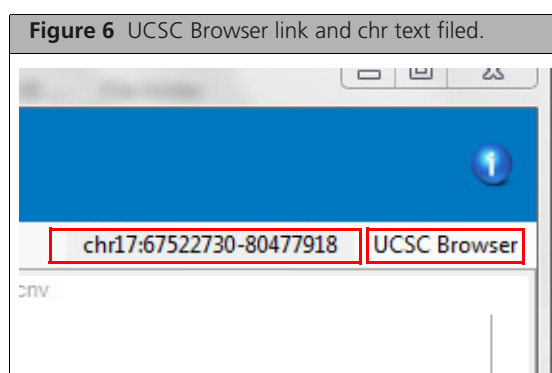
**To reset your customized view back to the default whole genome view:**

1. Click **Reset**.

## To use the CNV Viewer to investigate your copy number changes:

### Option #1

1. Use the *Zoom In* command or click, then drag your mouse cursor across a region of interest.
2. Once the Viewer has zoomed into a chromosome, a *UCSC Browser* button appears (upper right corner). (Figure 6)



3. Click on the **UCSC Browser** button.  
The UCSC website page appears (Figure 7) and displays the current region based on the chromosome positions listed in the *chr* text box. (Figure 6)

### Option #2

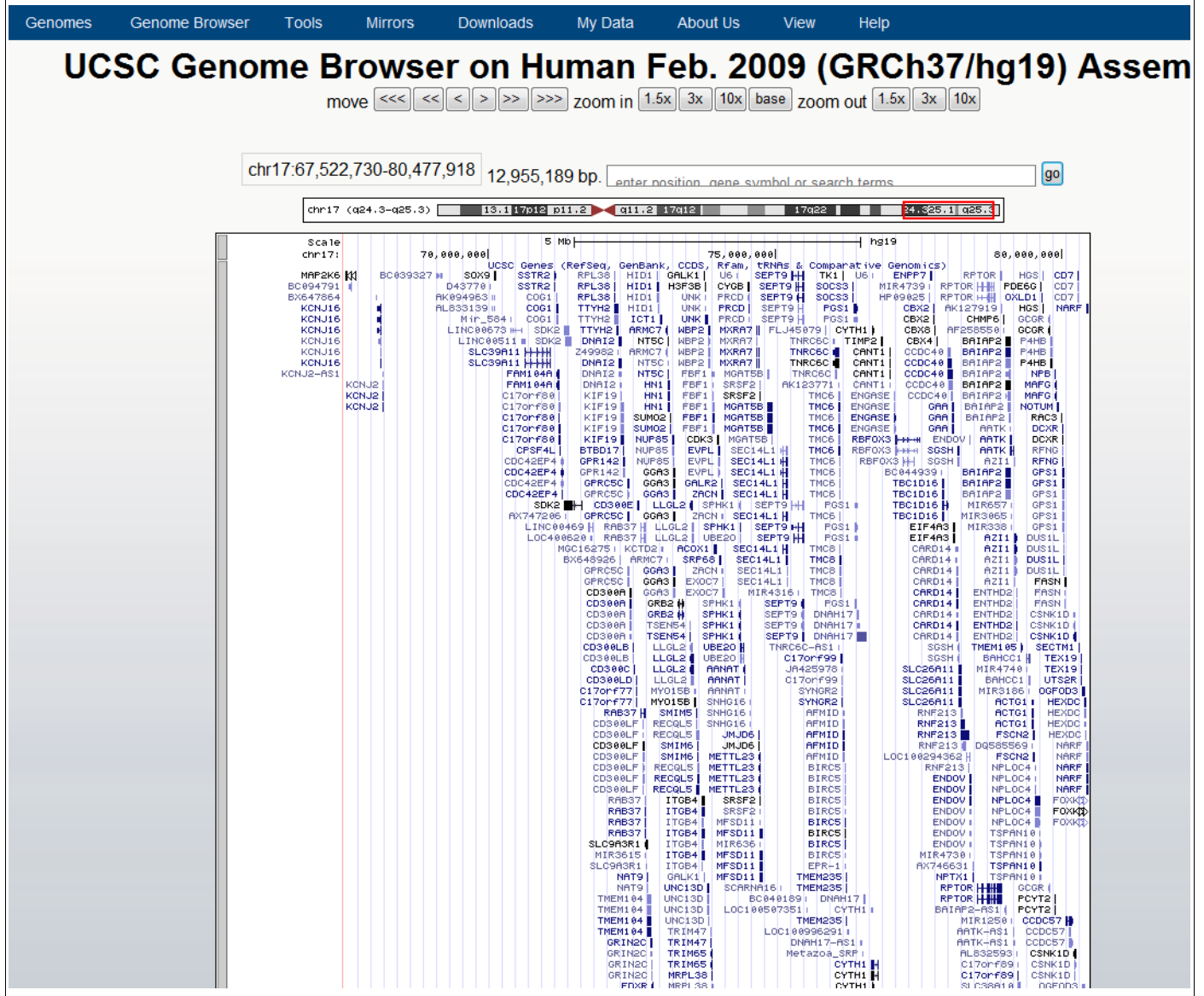
1. Click inside the *chr* text box (Figure 6), then manually enter your chromosome positions. You must use one of the following formats:
  - **chr17:67522730-80477918**
  - **chr17:67,522,730-80,477,918**
2. Press **Enter**.



**NOTE:** The region displayed in the CNV Viewer may be smaller than the chromosome positions you entered, because the CNV Viewer auto-adjusts your start and stop positions next to the nearest available start and stop markers.

3. (Optional) Click on the **UCSC Browser** button.  
The UCSC website page appears (Figure 7) and displays the current region based on the chromosome positions listed in the *chr* text box. (Figure 6)

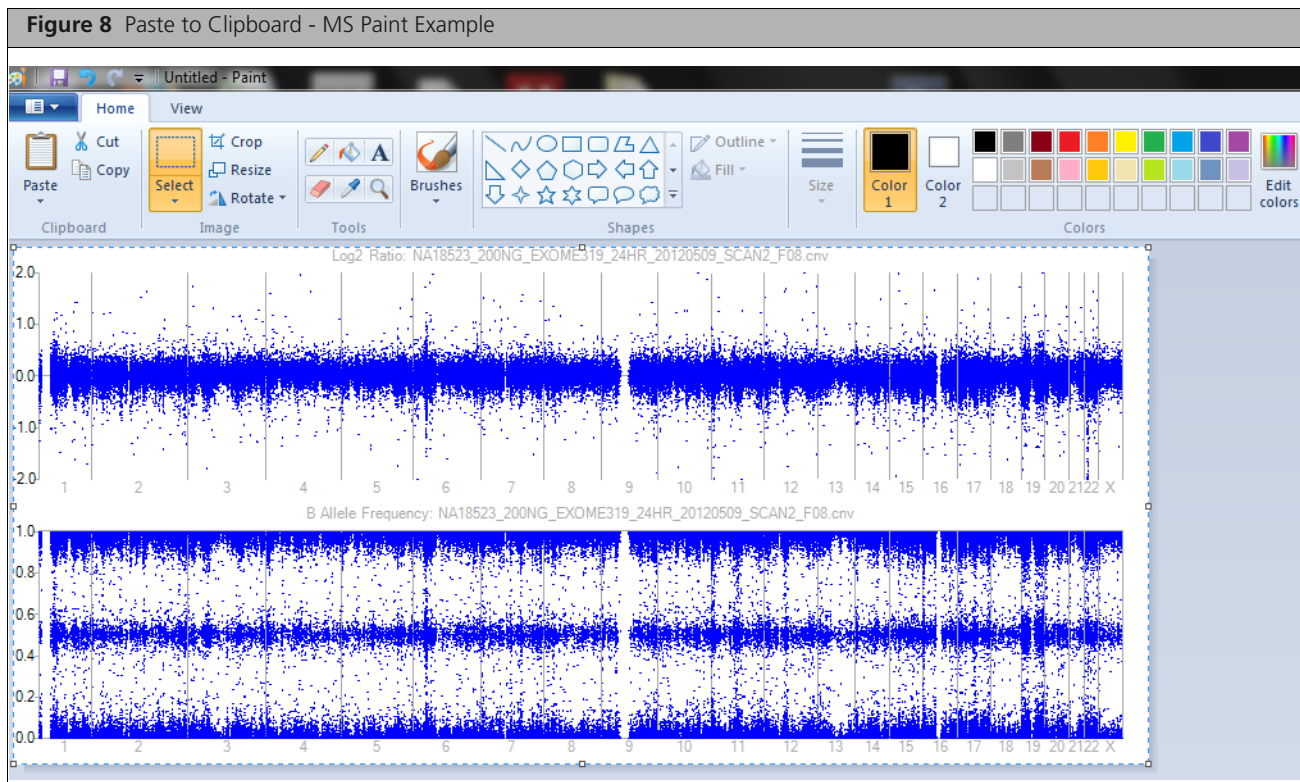
**Figure 7** UCSC website page





### To copy the current view to your Clipboard:

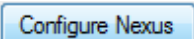
1. Click **Edit -> Copy to clipboard**.
2. Use the paste command (**Ctrl-V**) to copy the current view into another software application, such as MS Paint. (Figure 8)



## Further Copy Number Analysis Using BioDiscovery's Nexus Software

Use the newly generated **AxiomGT1.cnv.txt** and **<annot>.probemappings.txt** with BioDiscovery's Nexus software to perform copy number analysis.

**Do the following to configure Axiom CNV Summary Tool output data to work with BioDiscovery's Nexus software:**

1. Click the CNV Summary Tool's  button (bottom right).

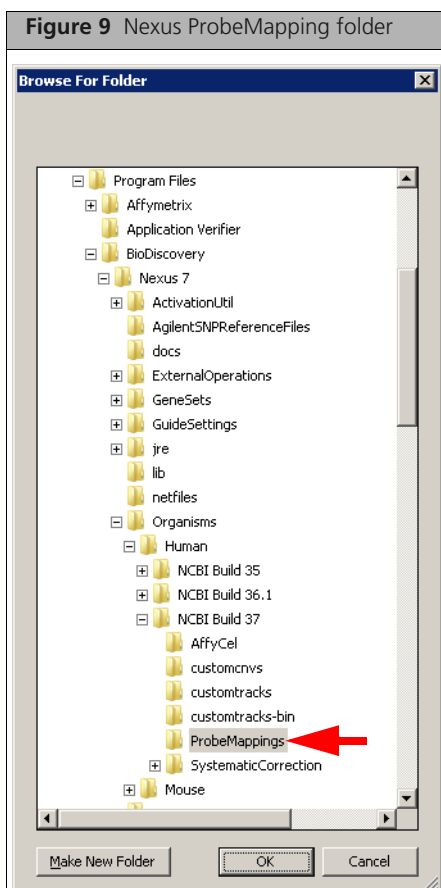


**NOTE:** If Nexus is not detected on your system, the *Configure Nexus* button is disabled.

If multiple versions of Nexus are detected, a drop-down menu appears. Use this menu to select the appropriate version of Nexus. This menu does not appear if only one version of Nexus is detected.

A file window appears.

2. Click to select the *probe mapping .txt* file. This file resides in your master Data Directory folder you setup earlier. See [Step 1 on page 8](#).
3. Click **Open**.  
An Explorer window appears. ([Figure 9](#))
4. Navigate Nexus's *ProbeMappings* directory/folder, then click **OK**.

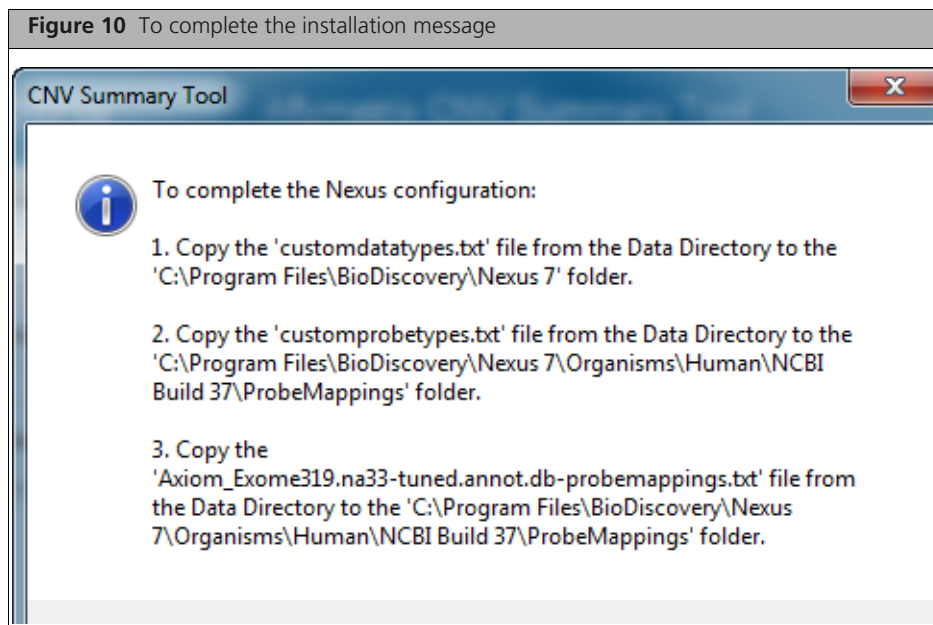


The message *Configuration Complete* appears.

5. Click **OK**.

**!** **IMPORTANT:** You are responsible for knowing the location of Nexus's *ProbeMappings* folder. If you are unsure of its location, contact BioDiscovery. In most cases, the Nexus Probe Mapping folder resides here: *C:\Program Files\BioDiscovery\NexusX*

If you do not have access (Administrator Privileges) to some of your computer's folders, a message with file copying instructions appears. (Figure 10) Follow the 3 steps shown to configure the Nexus software manually.



6. Use the Nexus software as you normally would.

If you want to perform a GC Correction in Nexus, see *Performing GC Correction in Nexus* on page 18

## Performing GC Correction in Nexus

**! IMPORTANT:** You must first download an appropriate BED file from [affymetrix.com](http://affymetrix.com).

**To download a BED file from [affymetrix.com](http://affymetrix.com):**

1. Go to [www.affymetrix.com](http://www.affymetrix.com).
2. Login as you normally would or click **Register**, then follow the on-screen instructions.
3. Click **Products** -> **Products** (top left).  
The Products page appears.
4. Click **Microarray Solutions** (left pane).  
The Microarray Solutions pane appears.
5. Under the **DNA Analysis Solutions** header, click to choose the option you want.  
Example: *Genome-Wide Genotyping for Human Disease Research*.  
For the *Genome-Wide Genotyping for Human Disease Research* example, 3 options appear.
6. Click the Arrays adjacent **[+]** button. (Figure 3.1)



For the *Genome-Wide Genotyping for Human Disease Research* example, the following page appears. (Figure 3.2)

Figure 3.2 Arrays menu

Home > Products > Microarray Solutions > DNA Analysis Solutions > **Genome-Wide Genotyping for Human Disease Research**

## Genome-Wide Genotyping for Human Disease Research

Affymetrix is a leading provider of microarray solutions for targeted and genome-wide applications. Our powerful portfolio includes arrays, reagents, instruments, and informatics tools that enable you to detect common and rare single nucleotide polymorphisms (SNPs), copy number variants, and other genetic variations that can contribute to complex diseases.

### Arrays

#### Cartridges

##### Human

- Genome-Wide Human SNP Array 6.0
- Human SNP Array 5.0, Made to Order
- Mapping 500K Array Set
- Mapping 100K Set, Made to Order
- Mapping 10K 2.0 Array, Made to Order

#### Plates

##### Human

- Axiom® Genotyping Solution
- Axiom® Exome Genotyping Arrays**
- Axiom® Biobank Genotyping Arrays
- Axiom® miRNA Target Site Genotyping Arrays
- Axiom® World Arrays
- Axiom® Genome-Wide Population-Optimized Human Arrays
- Axiom® Genome-Wide Human Origins 1 Arrays
- Axiom® myDesign™ Human Genotyping Arrays

Back to Top >

Reagents +

Instruments & Software +

- For this example, click the **Axiom® Exome Genotyping Arrays**. (Figure 3.2)  
The Axiom® Exome Genotyping Arrays page appears.
- Click on the **Technical Documentation** tab. (Figure 3.3)

Figure 3.3 Technical Documentation tab.

Product Description   **Technical Documentation**   Required/Related Products

- Scroll down and locate *NetAffx Alignment Files*, then (for this example) click on **Axiom Exome319 BED File**. (Figure 3.4)

Figure 3.4 Technical Documentation tab.

**NetAffx Alignment Files**

[Axiom Exome BED File \(4.3 MB, 3/28/12\)](#)

[Axiom Exome319 BED File \(3.2 MB, 3/6/13\)](#)

A Windows Explorer window appears.



**NOTE:** BED files for *Axiom myDesign* arrays are provided directly to you from Affymetrix (for each custom array designed).

---

10. Save the zip file to a convenient location.

**Do the following to submit your BED file to BioDiscovery for GC Correction:**

1. Extract the downloaded BED.zip file, then contact BioDiscovery and tell them you need a GC Correction file created from a BED file.

BioDiscovery will respond with an email containing a hyperlink.

2. Click on the hyperlink provided by BioDiscovery.

The following form appears. ([Figure 3.5](#))

3. Complete the form, click **Browse** to upload your unzipped BED file, then click **Submit File**.

Figure 3.5 BioDiscovery Form

### Create Systematic Correction File

This tool reads genome coordinates and creates a systematic correction file and sends email with the link to download the file

Enter email address:

Genome/Assembly:

Manufacturer:

Use Restriction Enzymes in fragmentation step:

Select neighborhood sizes to calculate GC%:

☐ 500b
 ☐ 1kb
 ☒ 4kb
 ☐ 5kb
 ☐ 10kb
 ☐ 50kb
 ☒ 100kb
 ☐ 150kb
 ☐ 500kb
 ☒ 1mb

Reject probes if fraction of unknown bases (N) exceeds:

Supported data formats.

- [Browser Extensible Data \(BED\)](#)
- Genomic Coordinate Position  
chrN:start-end

Paste in data:

Or upload data from a file:

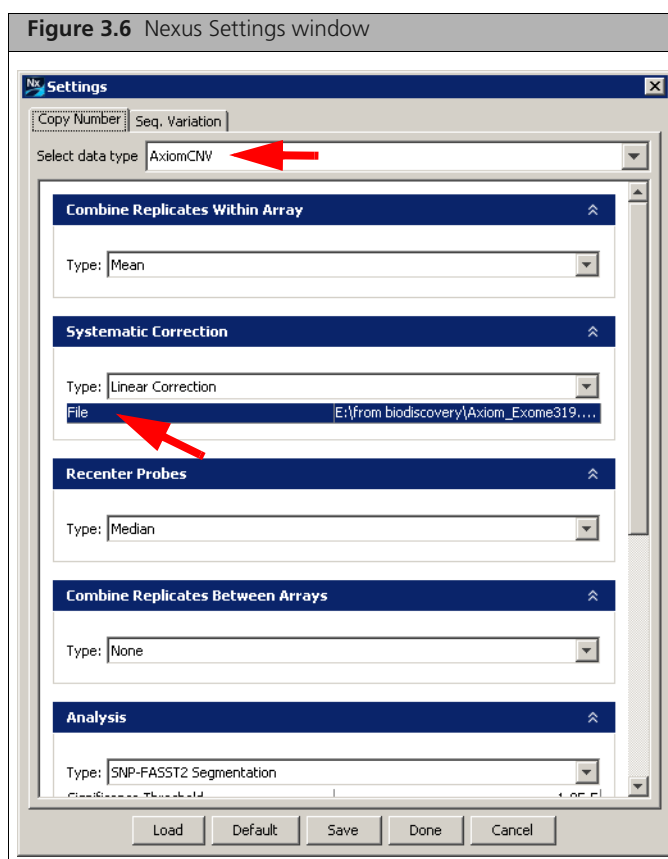
Axiom\_Exom....na33.bed



**NOTE:** BioDiscovery will email you a second hyperlink to download the GC Correction file for use with their Nexus software.

4. Click on the hyperlink to download/save the GC Corrected BED file. Make sure you save the file to a convenient location.
5. Open the *BioDiscovery Nexus* application as you normally would.
6. Click **Settings**.

The following window appears. (Figure 3.6)



7. From the *Select data type* drop-down menu, click **AxiomCNV**. (Figure 3.6)
8. From the *Systematic Correction* drop-down menu, select your GC Correction Type.
9. Click the *File* banner (Figure 3.6), then select your GC Correction file.
10. Use the other applicable drop-down menu selections to complete the Settings form, then click **Done**.
11. Use the Nexus software as you normally would.



## Calculations of log2 Ratios and B allele Frequencies

Affymetrix Axiom® Arrays, designed to detect genome-wide associations with SNPs and indels, can also detect copy number variations. The CNV detection method is based on computed log2 ratios and B allele frequencies (BAFs) at individual markers across the genome.

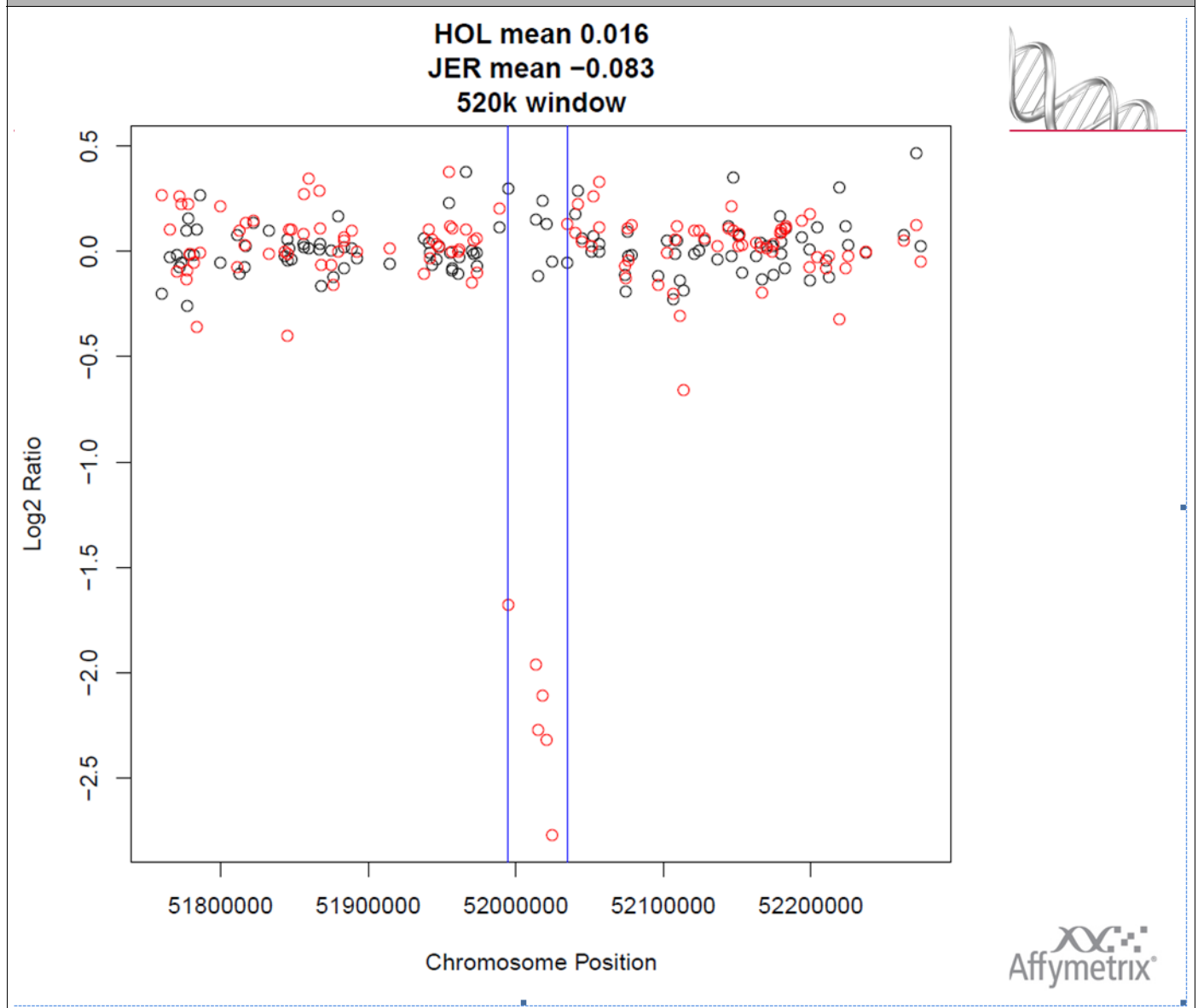
**The log2 ratios and B allele frequencies are calculated as follows:**

### Computing log2 Ratio

Log2 ratios are computed at each marker site as the sum of the A and B allele intensities for the marker normalized by the median intensity of that marker in phenotypically normal individuals, assumed to represent the normal copy number state at that marker site.

- a)** Total Intensity,  $T = \text{Intensity}_A + \text{Intensity}_B$
- b)** Across a reference set of samples for each marker determine a Reference value  $R = \text{median}(T_1, T_2, \dots, T_N)$ . If most individuals are expected have normal copy number states, one approach is to create the reference based on the  $N$  (~96) individuals genotyped on the plate. For the X chromosome, median values are taken over only Female samples. For the Y chromosome, median values are taken over only Male samples.
- c)** For each marker for each individual sample,  $\log_2 \text{ratio} = \log_2(T) - \log_2(R)$

The log2 ratio computed above may be ordered genomically and inspected visually for regions diverge from zero. [Figure 1](#) (below) shows a region with significant deletion in the genome of Jersey cattle relative to Holstein. The intensities were produced by genotyping with the Axiom® Genome-Wide BOS 1 Bovine array by Rincon et al [1]. ([Figure 1](#))

**Figure 1** log2 ratio values for copy number loss from Rincon, et al.

### B allele Frequencies (BAF)

The BAF at a marker is the ratio of the B allele intensity to the sum of the allele intensities. The raw BAFs are standardized based on mean BAFs for homozygous and heterozygous genotype calls for each marker and further scaled to be within [0,1]. (Figure 2)

**Figure 2** BAF marker equations

$$\text{Raw BAF } (\lambda) = \frac{\text{Intensity}_B}{\text{Intensity}_A + \text{Intensity}_B}$$

Standardized BAF =

$$\frac{\lambda - \lambda_{AB}}{|\lambda_{AA} - \lambda_{AB}|} \text{ when } \lambda < \lambda_{AB}$$

$$\frac{\lambda - \lambda_{AB}}{|\lambda_{BB} - \lambda_{AB}|} \text{ when } \lambda \geq \lambda_{AB}$$

where

$$\lambda_{AA} = \text{mean}(\lambda \mid AA)$$

$$\lambda_{AB} = \text{mean}(\lambda \mid AB)$$

$$\lambda_{BB} = \text{mean}(\lambda \mid BB)$$

For markers in the non-PAR regions of chromosome X, mean values are taken over only Female samples. For markers in the PAR regions of chromosome X, mean values are taken over all samples. For markers on the Y Chromosome, mean values are taken over only Male samples.

When there are no samples for a specific genotype at a marker site, then mean values must be interpolated.

If only  $\lambda_{AA}$  is missing, then  $\lambda_{AA} = \lambda_{AB} - (\lambda_{BB} - \lambda_{AB})$

If only  $\lambda_{AB}$  is missing, then  $\lambda_{AB} = (\lambda_{AA} + \lambda_{BB})/2$

If only  $\lambda_{BB}$  is missing, then  $\lambda_{BB} = \lambda_{AB} + (\lambda_{AB} - \lambda_{AA})$

If two mean values are missing, then the missing value is set to the median of the corresponding value of all probesets:

$$\lambda_{AA} = \text{median}(\lambda_{AA})$$

$$\lambda_{AB} = \text{median}(\lambda_{AB})$$

$$\lambda_{BB} = \text{median}(\lambda_{BB})$$

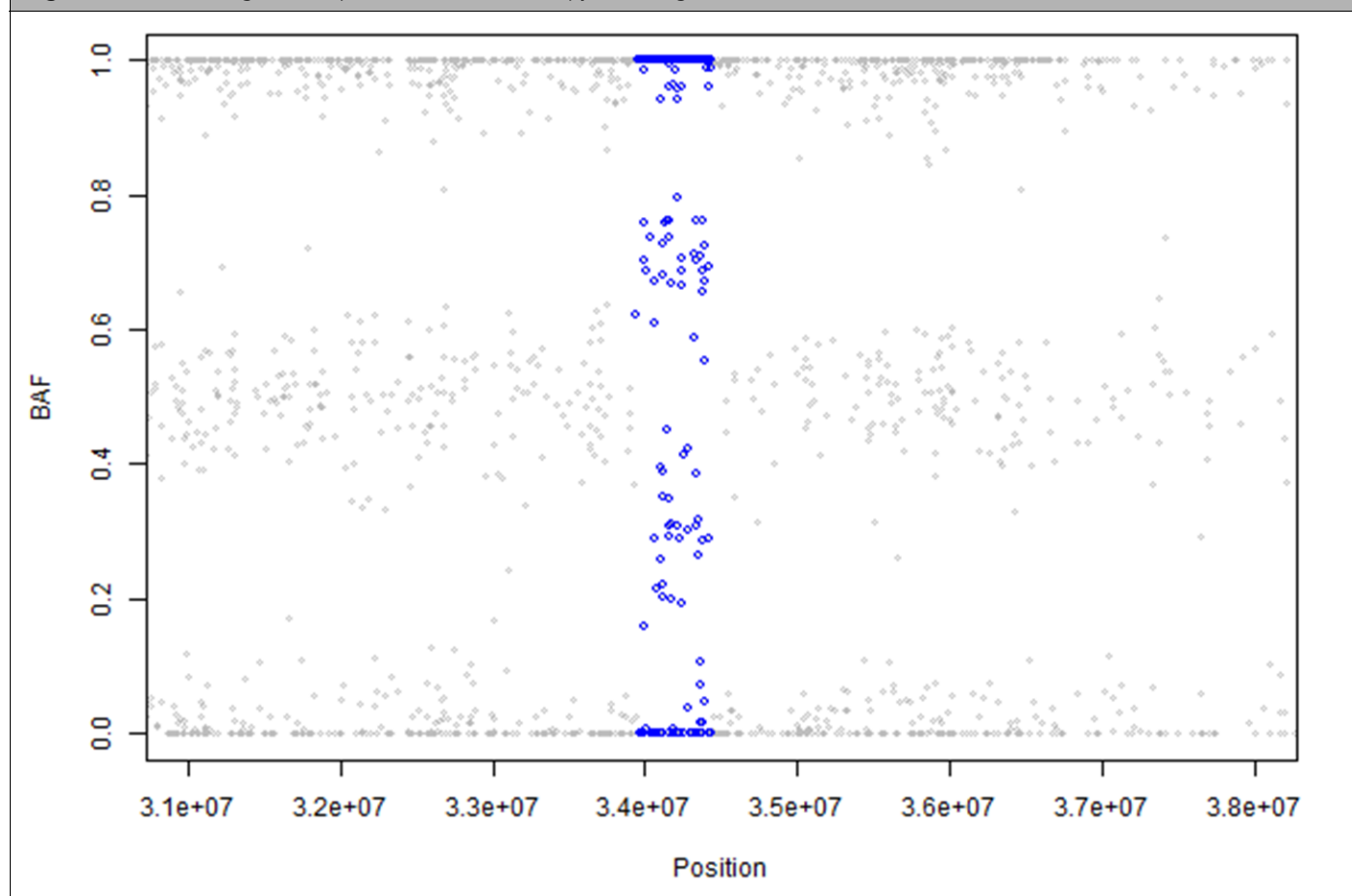
Scaled BAF =  $0.5 * (\text{Standardized BAF} + 1.0)$

The Scaled BAFs are truncated to be within [0,1]

The BAFs computed above may be ordered genomically and inspected visually for regions diverge from the expected pattern of 3 bands at 0, 0.5 and 1.0.

Figure 3 shows a region with a copy number 3 in the genome of a YRI HapMap sample. The intensities were produced by genotyping with Axiom® Genome Wide PanAFR array.

**Figure 3** BAFs showing a band split characteristic of a copy number gain (shown in blue).



## References

1. Rincon G, Weber K, Van Eenennaam A, et al. Hot topic: Performance of bovine high-density genotyping platforms in Holsteins and Jerseys. *J Dairy Sci.* 94:6116-6121 (2011).