

# Loss of Heterozygosity (LOH) Analysis Getting Started Guide





(LOH) Analysis

Loss of Heterozygosity

Getting Started Guide

**Getting Started** 

ï

Setting Up the Loss of Heterozygosity Analysis

2

Analyzing and Examining Results

3

Sorting Data and Evaluating Loss of Heterozygosity

4

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

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document. This document is believed to be complete and accurate at the time of publication. In no event shall Applied Biosystems be liable for incidental, special, multiple, or consequential damages in connection with or arising from the use of this document.

Notice to Purchaser: License Disclaimer.

Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Applied Biosystems, either expressly, or by estoppel.

GeneMapper Software has not undergone specific developmental validation for human identification applications. Human identification laboratories analyzing single-source or parentage samples which choose to use GeneMapper Software for data analysis should perform their own developmental validation studies.

#### TRADEMARKS:

Applied Biosystems, AB (Design), ABI PRISM, GeneMapper, LIZ and SNaPshot are registered trademarks and GeneScan and SNPlex are trademarks of Applied Biosystems or its affiliates in the U.S. and/or certain other countries.

This product includes software developed by the Apache Software Foundation (http://www.apache.org/). Copyright © 1999-2000 The Apache Software Foundation. All rights reserved.

This product includes software developed by the ExoLab Project (http://www.exolab.org/). Copyright 2000 © Intalio Inc. All rights reserved.

JNIRegistry is Copyright © 1997 Timothy Gerard Endres, ICE Engineering, Inc., http://www.trustice.com.

AFLP is a registered trademark of Keygene N.V.

Oracle is a registered trademark of Oracle Corporation.

All other trademarks are the sole property of their respective owners.

© Copyright 2009, Applied Biosystems. All rights reserved.

Part Number 4403621 Rev. A 04/2009

### **Contents**

	Preface	V
	How to Use This Guide	v
	How to Obtain More Information	vi
	How to Obtain Support	viii
Chapter 1	Getting Started	1
	About LOH Microsatellite Analyses	2
	About the Example Data	4
	LOH Microsatellite Analysis Workflow	6
	GeneMapper® Software Terms	7
	Starting the Software and Logging In	7
	Using This Guide With Your Own Sample Files	8
	Alternatives to the Procedures in This Guide	8
Chapter 2	Setting Up the Loss of Heterozygosity	
	Analysis	9
	Overview	10
	Creating a Kit, Panel, and Markers	11
	Creating a New Project and Adding Sample Files	15
	Setting Analysis Parameters and Table Settings for the Project	17
	Performing the Initial Analysis on the Project	
	Creating a Bin Set and Generating Bins	20
	Using the Auto Bin Feature	35

Chapter 3	Analyzing and Examining Results	45
	Editing the Analysis Method	46
	Analyzing the Project	48
	Examining the Results	48
Chapter 4	Sorting Data and Evaluating Loss of	
	Heterozygosity	53
	Overview	54
	Sorting the LOH Data	56
	Generating a Report to Calculate and Evaluate LOH	58
	Editing or Creating a Custom LOH Report Setting	59
Chapter 5	Printing and Exporting Results	65
	Printing Results	66
	Exporting Results	
	Index	69

### **Preface**

### How to Use This Guide

### Purpose of This Guide

The GeneMapper® Software Version 4.1 LOH Analysis Getting Started Guide provides brief, step-by-step instructions for sizing, genotyping, and evaluating LOH microsatellite data generated using any of the compatible Applied Biosystems electrophoresis instruments and Data Collection Software. It describes how to troubleshoot, print and export data, and create reports. It is designed to help you quickly learn to use basic functions of the GeneMapper Software.

Audience

This guide is intended for novice GeneMapper Software users.

### **Assumptions**

This guide assumes that:

- You have installed GeneMapper Software version 4.1 as described in the *GeneMapper® Software Version 4.1 Installation and Administration Guide* (PN 4403614).
- You have a working knowledge of the Microsoft® Windows® operating system.

### **Text Conventions**

This guide uses the following conventions:

- Bold indicates user action. For example:
   Type 0, then press Enter for each of the remaining fields.
- *Italic* text indicates new or important words and is also used for emphasis. For example:

Before analyzing, *always* prepare fresh matrix.

• A ▶ symbol separates successive commands you select from a drop-down or shortcut menu. For example:

Select File ▶ Open ▶ Spot Set.

Right-click the sample row, then select **View Filter > View All Runs**.

### User Attention Words

Two user attention words appear in Applied Biosystems user documentation. Each word implies a particular level of observation or action as described below:

**Note:** Provides information that may be of interest or help but is not critical to the use of the product.

**IMPORTANT!** Provides information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.

Examples of the user attention words appear below:

**Note:** The size of the column affects the run time.

**Note:** The Calibrate function is also available in the Control Console

**IMPORTANT!** To verify your client connection to the database, you need a valid Oracle user ID and password.

**IMPORTANT!** You must create a separate Sample Entry Spreadsheet for each 96-well plate.

### Safety Alert Words

Safety alert words also appear in user documentation. For more information, see the *GeneMapper® Software Version 4.1 Installation and Administration Guide* (PN 4403614).

### **How to Obtain More Information**

### Safety Information

For safety information, see the *GeneMapper® Software Version 4.1 Installation and Administration Guide* (PN 4403614).

### Software Warranty and License

For all warranty and licensing information, see the *GeneMapper*® *Software Version 4.1 Installation and Administration Guide* (PN 4403614).

### Related Documentation

The following related documents are shipped with the software:

- GeneMapper® Software Version 4.1 Installation and Administration Guide (PN 4403614) Provides procedures for installing, securing, and maintaining version 4.1 of the GeneMapper Software.
- GeneMapper® Software Version 4.1 Getting Started Guides for microsatellite analysis (PN 4403672), loss of hetereozygosity (LOH) analysis (PN 4403621), AFLP® system analysis (PN 4403620), SNaPshot® kit analysis (PN 4403618), and SNPlex™ system analysis (PN 4403617) − Five guides that explain how to analyze the application-specific example data provided with the GeneMapper Software. The guides provide brief, step-by-step procedures for the analysis of microsatellite, LOH, AFLP® system, SNaPshot® kit, and SNPlex™ system data generated by compatible Applied Biosystems electrophoresis instruments and Data Collection Software. The guides are designed to help you quickly learn to use basic functions of the GeneMapper Software.
- GeneMapper® Software Version 4.1 Online Help Describes the GeneMapper Software and provides procedures for common tasks. Access online help by pressing F1, selecting Help
   Contents and Index, or clicking in the toolbar of the GeneMapper window.
- GeneMapper® Software Version 4.1 Quick Reference Guide
   (PN 4403615) Provides workflows for specific analysis types
   and lists instruments, software, and analysis applications
   compatible with the GeneMapper Software.
- GeneMapper® Software Version 4.1 Reference and Troubleshooting Guide (PN 4403673) Provides reference information such as theory of operation and includes troubleshooting information.

Portable document format (PDF) versions of this guide and the other documents listed above are available on the *GeneMapper Software Version 4.1 Documentation CD*.

**Note:** For additional documentation, see "How to Obtain Support" on page viii.

### Obtaining Information from Online Help

The GeneMapper Software features an online help system that describes how to use each feature of the user interface. Access online help by pressing F1, selecting Help > Contents and Index, or clicking in the toolbar of the GeneMapper window.

### Send Us Your Comments

Applied Biosystems welcomes your comments and suggestions for improving its user documents. You can e-mail your comments to:

techpubs@appliedbiosystems.com

### **How to Obtain Support**

For the latest services and support information for all locations, go to <a href="http://www.appliedbiosystems.com">http://www.appliedbiosystems.com</a>, then click the link for Support.

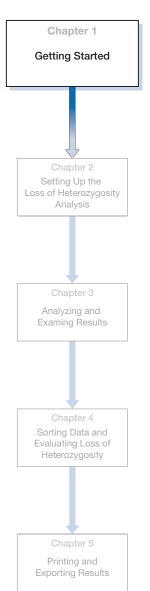
At the Support page, you can:

- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support
- Order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents
- Download PDF documents
- Obtain information about customer training
- Download software updates and patches

In addition, the Support page provides access to worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.



### **Getting Started**



This chapter includes:	
■ About LOH Microsatellite Analyses	2
About the Example Data	4
■ LOH Microsatellite Analysis Workflow	6
■ GeneMapper® Software Terms	7
■ Starting the Software and Logging In	7
■ Using This Guide With Your Own Sample Files	8
■ Alternatives to the Procedures in This Guide	8

### **About LOH Microsatellite Analyses**

### Microsatellite Markers

Microsatellite markers, also known as short tandem repeats (STRs), are polymorphic DNA loci consisting of a repeated nucleotide sequence. The repeat sequence can be from 2 to 7 base pairs long. The number of repeat units varies in a population, thereby creating multiple alleles for a microsatellite locus.

### Microsatellite Analysis

In a typical microsatellite analysis, microsatellite loci are amplified by PCR using fluorescently labeled forward and unlabeled reverse primers. The PCR amplicons are separated by size using electrophoresis; then the dye labeled products are identified by fluorescence detection. You can then use the GeneMapper Software to size and genotype the alleles.

#### What is LOH?

In the two-hit model used to describe inactivation of tumor suppressor genes (TSGs), the first mutation or "hit" results in a heterozygous state for the TSG with one wild-type and one mutant allele. If this is followed by a second "hit," in which all or part of the chromosome that contains the wild-type allele of the TSG is deleted, the chances of tumorigenesis increase. This phenomenon is known as Loss of Heterozygosity or LOH. While this is a rare event, it occurs more frequently in familial forms of cancer, in which a mutation of one of the TSG alleles is inherited.

### LOH Microsatellite Analysis

An LOH microsatellite analysis is the screening of tumor samples for LOH using microsatellite markers. Because LOH can be caused by deletion of genomic DNA regions containing the wild-type copy of a TSG, researchers can use microsatellite markers to screen samples for LOH.

A typical LOH assay compares amplified microsatellite markers from the suspected cancerous tissue to the same markers from the healthy tissue, both from the same individual. The healthy tissue should show two alleles for a given heterozygous microsatellite marker. If a given tumor sample has undergone LOH, it will be evident by a decrease in peak height of one of the two alleles (relative to the healthy allele peak heights) (Figure 1-1). The reason the tumor sample shows a decrease in peak height (instead of an absence of the peak) is that during isolation, the tumor sample is contaminated with healthy cells, thus introducing some wild-type DNA into the tumor specimen.

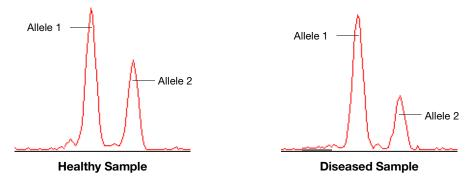


Figure 1-1 Diseased sample shows loss in peak height of Allele 2

After performing a microsatellite analysis, you can use the Report Manager in the GeneMapper® Software to calculate and compare peak height ratios of microsatellite alleles between healthy and diseased tissues, as well as identify and flag LOH candidates based on set peak-height ratio thresholds. You can alter these thresholds based on the observed level of wild-type DNA contamination in the tumor samples.

#### **Custom Primers**

Applied Biosystems provides custom primers for PCR amplification of LOH microsatellite markers. For more information, visit the Applied Biosystems Web site at www.appliedbiosystems.com.

### **Compatible Instruments**

For information about Applied Biosystems electrophoresis instruments that are compatible with LOH microsatellite analyses, see the *GeneMapper*<sup>®</sup> *Software Version 4.1 Quick Reference Guide* (PN 4403615).

### **About the Example Data**

### Sample File Naming Conventions

The naming convention for the sample files used in this guide allows you to take advantage of the LOH Default report setting provided with the GeneMapper® Software. Specifically, healthy and tumor samples from the same individual start with the same letter or number, so they list consecutively when sorted by sample file and marker.

**Example:** 1\_Healthy.fsa, 1\_Tumor.fsa, 2\_Healthy.fsa, 2\_Tumor.fsa, and so on

When analyzing your own LOH data, if you want to use the LOH Default report setting, your sample files must be named as described above.

If your LOH data sample files are not named as described above, you can still use the sorting and LOH reporting features in the GeneMapper® Software, but you will have to edit the report setting as described in "Editing or Creating a Custom LOH Report Setting" on page 59. Below is an example naming convention in which healthy and tumor samples from the same individual do *not* list consecutively after being sorted by sample file and marker.

**Example:** Healthy\_1.fsa, Healthy\_2.fsa, ...,Tumor\_2.fsa, Tumor\_1.fsa, and so on

#### Sample File Location

To perform the exercise described in this getting started guide, use the four sample files (.fsa) located on your computer hard drive at:

<drive>:\AppliedBiosystems\GeneMapper\Example
Data\LOH

**Note:** The above location will vary depending on the installation of the GeneMapper<sup>®</sup> Software. The default installation is the D drive.

### Instrument and Size Standard

Sample files were generated by running PCR-amplified and fluorescently tagged LOH samples on an ABI PRISM® 3100 Genetic Analyzer using the GeneScan<sup>TM</sup> 500 LIZ® Size Standard.

### Marker Information

The sample files contain the following six markers.

Marker	Dye Label	Allele Size Range (bp)
R5	Blue	120 – 186
R26	Blue	193 – 295
R14	Blue	300 – 470
R21	Yellow	100 – 160
R24	Yellow	170 – 250
R2	Green	157 – 204

You will use this marker information when creating a panel and markers in Chapter 2, "Setting Up the Loss of Heterozygosity Analysis."

### **LOH Microsatellite Analysis Workflow**

The following flowchart summarizes the steps for performing an LOH analysis using the GeneMapper<sup>®</sup> Software:

### Set Up the Loss of Heterozygosity Analysis (Chapter 2)

- 1. Create a kit, panel, and markers for the project.
- 2. Create a new project and add sample files.
- 3. Set the analysis parameters and table settings for the project.
- 4. Perform an initial analysis.
- Create a bin set and generate bins (using Auto Bin).

### Analyze and Examine Results (Chapter 3)

- 1. Edit the analysis method to specify a bin set.
- 2. Analyze the samples in the project.
- 3. Examine the results.

### Sort Data and Evaluate Loss of Heterozygosity (Chapter 4)

- 1. Sort the LOH data in the Genotypes tab by sample file and marker.
- 2. Generate a report to calculate and evaluate LOH.
- 3. (Optional) Edit or create a custom LOH report setting.

### Print and Export Results (Optional) (Chapter 5)

- · Print results.
- · Export results.

### GeneMapper® Software Terms

Term	Definition
analysis parameters	A collection of user-defined settings (including an analysis method, size standard, and panel) that determine the sizing and genotyping algorithms used by the GeneMapper <sup>®</sup> Software to analyze all sample files in a project.
bin	A fragment size (bp) and dye color that define an allele within a marker. You create a bin for each possible allele associated with a marker.
bin set	A collection of bins (allele definitions), typically specific to a set of experimental conditions.
marker	A microsatellite marker is defined by a name, fragment size range (bp), dye color, and repeat length.
panel	A group of markers. In the GeneMapper Software, you associate a panel with a bin set to provide bin definitions for the markers.
kit	A group of panels.

### Starting the Software and Logging In

To start the GeneMapper® Software and log in:

- 1. Select Start ▶ All Programs ▶ Applied Biosystems ▶ GeneMapper ▶ GeneMapper 4.1.
- **2.** In the Login to GeneMapper dialog box:
  - **a.** Type the **User Name** and **Password** assigned by your system administrator.
  - b. Click OK.

### Using This Guide With Your Own Sample Files

In addition to using this guide to analyze the example data provided with the software, you can use this guide to lead you through the general LOH microsatellite analysis workflow when analyzing your own sample files. For information on advanced software features, see the *GeneMapper® Software Online Help*.

### Alternatives to the Procedures in This Guide

#### Overview

This guide presents one of several possible solutions for analyzing microsatellite LOH data using the GeneMapper<sup>®</sup> Software. Once you have completed the exercises in this document, you will most likely want to tailor the process to fit the requirements of your laboratory. This section provides you with a summary of several alternatives and where to go for further information.

### Using Autoanalysis to Set Up Projects

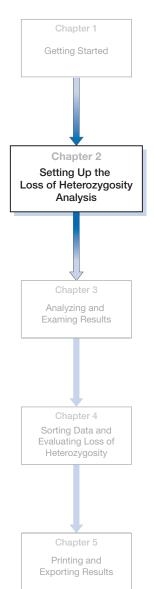
The GeneMapper Software includes an Autoanalysis feature that can eliminate most of the tasks leading up to the analysis of a microsatellite LOH project. Much of Chapter 2, "Setting Up the Loss of Heterozygosity Analysis," explains how to manually create, add samples to, and analyze projects for use in microsatellite LOH projects. When configured for Autoanalysis, the GeneMapper Software can automatically accomplish these tasks by coordinating with the Data Collection Software. For a more detailed explanation of how to use the Autoanalysis feature to set up microsatellite LOH projects, see the *GeneMapper® Software Version 4.1 Installation and Administration Guide* (PN 4403614).

### Using the Command Line Interface to Set Up Projects

The GeneMapper Software features a command line interface that can perform most of the major functions of the software. The command line interface can be a useful tool when analyzing microsatellite LOH projects because it automate many of the tasks explained in Chapter 2, "Setting Up the Loss of Heterozygosity Analysis." For a complete description of the command line interface and how it can be used to automate the functions of the GeneMapper Software, see the GeneMapper® Software Version 4.1 Installation and Administration Guide (PN 4403614).

2

## Setting Up the Loss of Heterozygosity Analysis



This chapter includes:	
Overview	10
■ Creating a Kit, Panel, and Markers	11
■ Creating a New Project and Adding Sample Files	15
■ Setting Analysis Parameters and Table Settings for the Project	17
Performing the Initial Analysis on the Project	26
■ Creating a Bin Set and Generating Bins Using the Auto Bin Feature	35

### **Overview**

### In This Chapter

In this chapter you will learn how to:

- Create a kit, panel, and markers
- Create a new project and add sample files
- Set analysis parameters and display settings for the project
- Perform the initial analysis on the project
- Create a bin set and generate bins using the Auto Bin feature

### For More Information

This chapter contains basic procedures. It does not describe all features and parameters in the GeneMapper Software. For more detailed information on topics presented in this chapter, see the following topics in the *GeneMapper*® *Software Online Help*:

- Creating a New Kit
- Creating a Custom Panel
- · Creating Markers
- · Creating a Project
- · Adding Samples
- Applying Analysis Settings
- Starting Analysis
- Creating a New Bin Set
- Using the Auto Bin Function

Online help is available from the Help menu, by clicking  $\bigcirc$ , or by pressing  $\mathbf{F1}$ .

### Creating a Kit, Panel, and Markers

#### Overview

You create the following hierarchical objects in the Panel Manager:

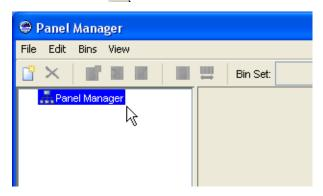
- **Kit** A group of panels
- **Panel** A group of markers
- Marker A fragment size range (bp), dye color, and repeat length

**Note:** In this guide, you will learn how to create panels and markers. However, you can also import panels (text files) that contain marker information. For example, panel files are available in the GeneMapper<sup>®</sup> Software for some of the LMS kits available from Applied Biosystems. For more information on importing panels, see the *GeneMapper*<sup>®</sup> *Software Online Help*.

### Creating a Kit, Panel, and Markers

To create a kit, panel, and markers:

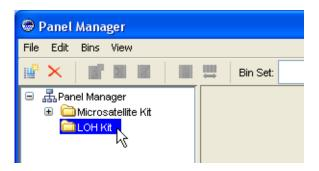
- 1. Open the Panel Manager by clicking ☐ (Tools ▶ Panel Manager).



**3.** In the New Kit dialog box, type **LOH Kit** for the Kit Name, select **Microsatellite** for the Kit Type, then click **OK**.

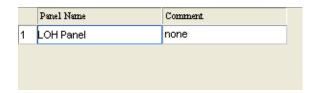


LOH Kit appears in the Navigation Pane (left side).



4. Select the LOH Kit in the Navigation Pane, then click (File ▶ New Panel).

**5.** In the right pane of the Panel Manager, select **New Panel**, type **LOH Panel** for the Panel Name, then press **Enter**.



LOH Panel appears under LOH Kit in the Navigation Pane (left side).



- 6. Select the LOH Panel in the Navigation Pane, then click (File ▶ New Marker).
- **7.** In the right pane of the Panel Manager, type the following marker information:

	Marker Name	Dye Color	Min Size	Max Size	Control Alleles	Marker Repeat
1	R5	blue	120.0	186.0		2

**8.** Repeat steps 6 through 7 to create the following markers:

Marker	Dye Color	Minimum Size (bp)	Maximum Size (bp)	Marker Repeat
R26	Blue	193	295	2
R14	Blue	300	470	2
R21	Yellow	100	160	2
R24	Yellow	170	250	2
R2	Green	157	204	2

**9.** Click **OK** to apply your changes and close the Panel Manager.

### **Next Steps**

Create a new project and add sample files (.fsa) to it as described on page 15.

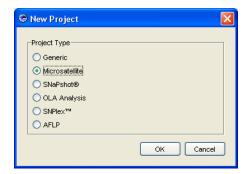
### Creating a New Project and Adding Sample Files

#### Overview

You create a project and add samples to the project in the GeneMapper window.

Creating a New Project and Adding Sample Files To create a new project and add sample files:

1. Click (File ▶ New Project).



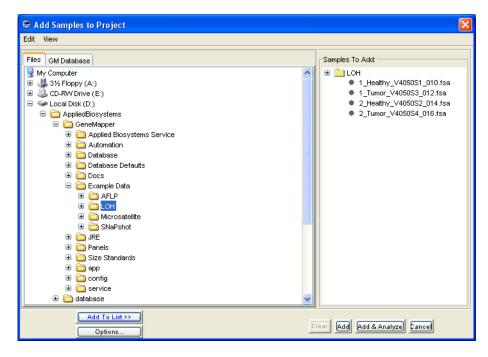
- 2. In the New Project dialog box, select **Microsatellite**, then click **OK**
- **4.** In the Add Samples to Project dialog box, in the Files tab, navigate to:

**Note:** The above location will vary depending on the installation of the GeneMapper<sup>®</sup> Software. The default installation is the D drive.

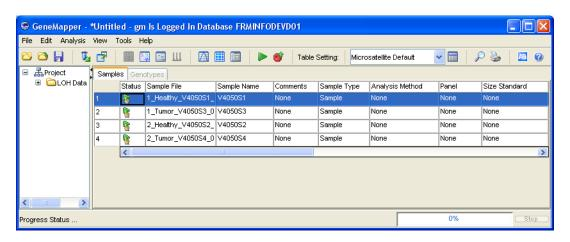
5. Select the LOH folder, click Add to List, then click Add.

**Note:** For this guide you added all sample files in the LOH folder. However, you can add a subset of files from a folder by expanding the folder in the left pane, pressing and holding Ctrl, then selecting individual files before clicking Add To List.





The four sample files from the LOH Data folder appear in the Samples tab, along with information entered in the Data Collection Software on the compatible Applied Biosystems electrophoresis instrument.



**Next Steps** Set analysis parameters and display settings for the project as described on page 17.

### Setting Analysis Parameters and Table Settings for the Project

#### Overview

You set analysis parameters and display settings for the project in the GeneMapper window.

Analysis parameters include:

- Analysis method (including bin set)
- Panel (set of markers)
- · Size standard

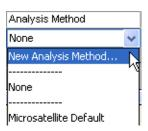
You set analysis parameters that determine the peak detection, sizing, and genotyping algorithms used by the GeneMapper® Software to analyze all sample files in a project.

Display settings include Table Settings and Plot Settings.

### Setting Analysis Parameters

To set analysis parameters for the project:

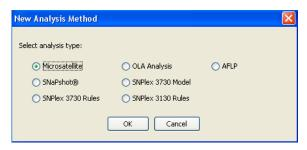
- **1.** Select the **Samples** tab in the GeneMapper window.
- 2. Click the first row in the Analysis Method column, then select New Analysis Method from the drop-down list.



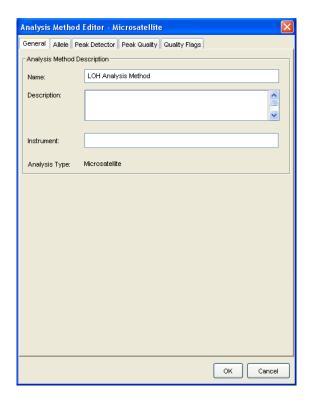
**Note:** You can also create a new analysis method from the Analysis Method tab in the GeneMapper Manager.



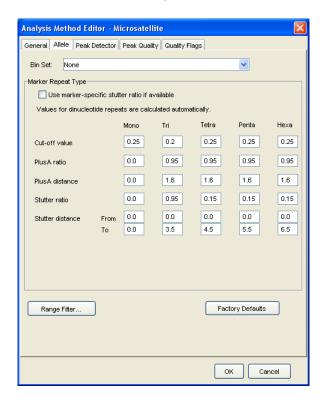
3. In the New Analysis Method dialog box, select Microsatellite for Analysis Type, then click OK.



- **4.** In the Analysis Method Editor dialog box, select and edit the five tabs:
  - General This tab includes reference information about the method. Type **LOH Analysis Method** for the Name. Optionally, type a description and the instrument on which the data was generated.

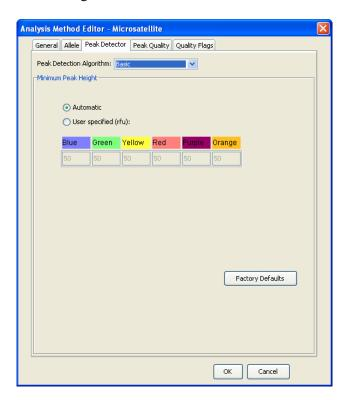


• Allele – This tab includes settings that determine allele calling. Select **None** for the Bin Set. Leave the default values for all other settings.





• **Peak Detector** – This tab includes settings that determine peak detection and sizing of peaks. Select **Basic** for the Peak Detection Algorithm. Leave the default values for all other settings.



• Peak Quality – This tab includes settings that determine when specific PQVs are left green (Pass) or flagged yellow (Check).

Type **2.0**. for the Max peak width (basepairs). Leave the default values for all other settings.





#### • Quality Flags – This tab includes:

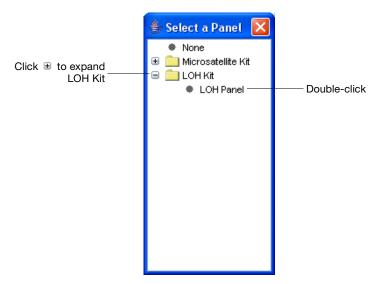
- Settings that determine the importance of individual flagged Process Quality Values (PQVs) to the overall Genotype Quality (GQ). You can weight each PQV from 0 to 1, with 0 being of no importance and 1 meaning very important.
- Threshold settings that determine when the SQ and GQ are flagged as Pass , Check , or Low Quality .
   The SQ and GQ are given initial scores of 1. The value of any flagged PQVs are then subtracted from 1 to give the final SQ and GQ scores.
- Assume Linearity Range, where the size calling algorithm assumes the fragment migration is linear for a given size range when calculating the Sizing Quality (SQ).

Leave the default values for all settings.



For details on analysis method parameters, see the *GeneMapper*® *Software Online Help*.

- **5.** Click **OK** to save the method and close the Analysis Method Editor dialog box.
- 6. Select the first row in the Panel column. From the Select a Panel dialog box, expand the LOH Kit, then double-click LOH Panel. (This is the Panel you created on page 11.)



7. Select the first row in the Size Standard column, then select GS500(-35, -250, -340)LIZ from the drop-down list.

**Note:** In the GeneMapper software, the following size standards are available for use with samples run with the GeneScan<sup>TM</sup> 500 LIZ<sup>®</sup> Size Standard:

- GS500(-250)LIZ excludes the 250-bp peak
- GS500(-35,-250,-340)LIZ excludes the 35-, 250-, and 340-bp peaks
- GS500LIZ includes all peaks present in the actual GeneScan™ 500 LIZ® Size Standard

Depending on your instrument, polymer type, and primer, it may be appropriate to choose one of the other size standards that omits peaks. Specifically, the 35-bp peak can be eclipsed by the neighboring primer peak, or the 250- and 340-bp peaks can migrate abnormally on the capillary electrophoresis instrument. Additionally, you can create your own custom size standards. For information on creating custom size standards, see the *GeneMapper*® *Software Online Help*.

- **8.** Fill down your selections to all sample rows in the Samples tab:
  - **a.** Click-drag across the Analysis Method, Panel, and Size Standard column headers to highlight all rows in all three columns.

Analysis Method	Panel	Size Standard
LOH Analysis Method	LOH Panel	GS500(-35,-250,-340)LIZ
None	None	None
None	None	None
None	None	None

**b.** Select **Edit** ▶ **Fill Down** (or press **Ctrl-D**).

### Selecting Table Setting

At the top of the GeneMapper window, select **Microsatellite Default** from the Table Settings drop-down list.

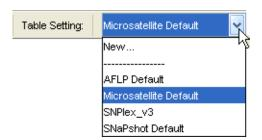


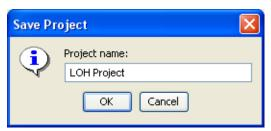
Table Settings control the information displayed in the Samples tab and Genotypes tab after analysis. Microsatellite Default is one of the default Table Settings provided with the GeneMapper Software.

You can also edit and create custom Table Settings in the GeneMapper Manager. For more information, see the *GeneMapper*® *Software Online Help*.

### Saving the Project

### To save the project:

- **2.** In the Save Project dialog box, type **LOH Project**, then click **OK**.



LOH Project appears in the title bar of the GeneMapper window.

**Next Steps** Perform the initial analysis on the project as described on page 26.

### Performing the Initial Analysis on the Project

#### Overview

Now that you have added sample files to and set analysis parameters for the project, perform an initial analysis to size the data so you have sample files available as reference data to create bins (allele definitions).

**Note:** Because you selected a Panel for the sample files in the project, the GeneMapper<sup>®</sup> Software will not only size the data but also try to genotype the data. However, because you did not specify a bin set in the analysis method, the Genotype Quality (GQ) will fail.

To perform the initial analysis:

- Analyze the project
- Review the SQ and contributing PQVs
- Examine the size standard
- View sample information (including raw data)
- Viewing samples plots

### Analyzing the Project

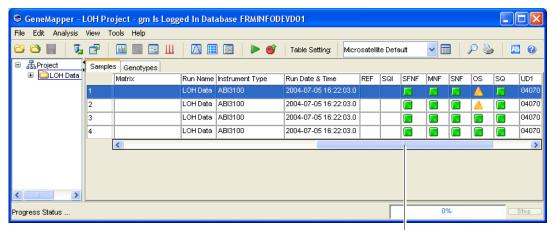
Click ► (Analysis ➤ Analyze).

The GeneMapper Software analyzes each sample in the project, displaying its progress in the Status Bar (lower left) of the GeneMapper window.

# Reviewing the SQ and PQVs

#### To review the Size Quality (SQ) and contributing PQVs:

- **1.** Make sure "Analysis Completed" appears in the Status Bar (lower left) of the GeneMapper window.
- **2.** Review the SQ by scrolling to the right in the Samples tab.



Click-drag scroll bar to right to view SQ column

If you followed the procedures and used the example data indicated in this guide, the SQ for each sample is (Pass). Most of the Process Quality Values (PQVs) that contribute to the SQ (SFNF, MNF, SNF, and OS) should also be .

Investigating Yellow A and Red SQs

IMPORTANT! When analyzing your own data, you may find the SQ to be ▲ (Check) or ♠ (Low Quality) and associated PQVs (SFNF, MNF, SNF, and OS) to be ▲, indicating issues with the size standard, data, or analysis parameters. To investigate and correct these issues, see "Examining the Size Standard" on page 28.

Note: Click of to sort the samples by SQ score. Samples with a SQ will be listed at the top of the Samples tab.

# Examining the Size Standard

#### To examine the size standard:

- Select all samples in the Samples tab by selecting Edit ➤ Select All.

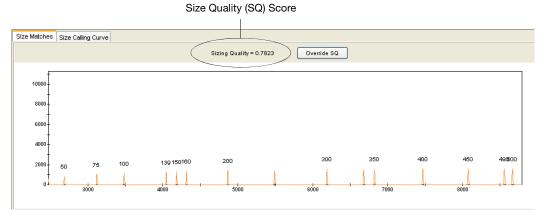


Figure 2-1 Size Match Editor - Size Matches tab

- **3.** Click the **Size Matches** tab to view the following for the selected sample:
  - Size Quality (SQ) score
  - · Size standard peaks
  - Size standard peak labels
- 4. Note the Sizing Quality score (Figure 2-1) for the sample. This score reflects how well the data from the size standard match the size standard you selected in the software. This score determines whether the SQ displays (Pass), (Check) or (Low Quality).

If you followed the instructions in this guide, the Sizing Quality is > 0.75 and the SQ displays (Pass).

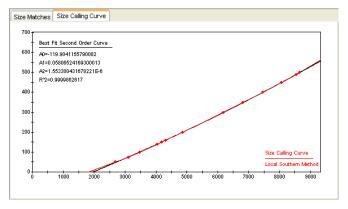
However, when analyzing you own data you may notice the Sizing Quality is less and the SQ displays △ (Check) or (Low Quality). For troubleshooting help, see Table 2-1 on page 30.

**5.** Determine if all peaks in the size standard are present and labeled correctly.

If you followed instructions in this guide, all peaks are present and labeled correctly as shown in Figure 2-1.

However, when analyzing you own data you may find some size standards peaks to be incorrectly labeled or missing. For troubleshooting help, see Table 2-1 on page 30.

**6.** Click the **Size Calling Curve** tab to view the size standard curve for the selected sample. You will see red data points representing the fragments from the size standard and a black best-fit curve.



- **7.** Select another sample from the left pane of the Size Match Editor, then repeat steps 3 through 6.
- **8.** Click **OK** to close the Size Match Editor.

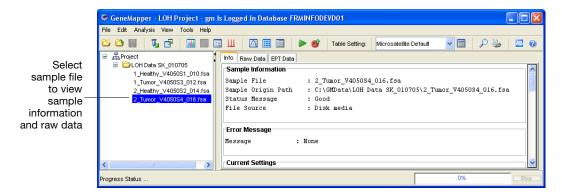
Table 2-1 Troubleshooting the size standard

Problem	Action				
Sizing Quality score is low and the SQ displays (Check) or (Low Quality), but all size standard peaks are present and labeled correctly.	Override the Sizing Quality by clicking <b>Override SQ</b> at the top of the Size Matches tab (Figure 2-1). Overriding changes the Sizing Quality score to 1.0, indicating the user verified the size standard.				
Some size standard peaks are not labeled correctly.	Edit, delete, and add size labels in the Size Matches tab, then click <b>Apply</b> to reanalyze the data with the updated sizing information. For more information, see the <i>GeneMapper® Software Online Help</i> .				
Some size standard peaks are not present.	Create a custom size standard in the software. For more information, see the GeneMapper® Software Online Help.				

For additional help in troubleshooting sizing problems, see the GeneMapper® Software Reference and Troubleshooting Guide.

#### Viewing Sample Information

To view information and raw data associated with individual sample files, select a sample file in the Navigation Pane (left), then select the Info or Raw Data tabs.

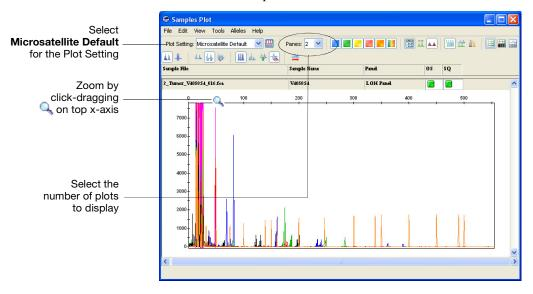


#### Viewing Sample Plots

To view the plots of the samples:

- **1.** Select **View > Samples** to display the Samples tab.
- 2. Select a sample (row) in the Samples tab. To select multiple samples, press and hold **Shift** or **Ctrl**. To select all samples, select **Edit** > **Select All**.
- 3. Click (Analysis ➤ Display Plots).

The Samples Plot window displays an electropherogram for each selected sample.



**4.** Select **Microsatellite Default** for the Plot Setting.

**Note:** Plot Settings control the information displayed in the Samples Plot window after analysis. Microsatellite Default is one of the default Plot Settings provided with the GeneMapper Software. You can also edit and create custom Plot Settings in the GeneMapper Manager. For more information, see the *GeneMapper® Software Online Help*.

#### **5.** Zoom on the x- and y-axes in the Samples Plot:

То	Then
Zoom on a specific region of the x-axis	Place the cursor on the top x-axis, then click-drag the q right or left to zoom all plots. Press and hold <b>Shift</b> while click-dragging to zoom only the selected plot.
	or
	Right-click the top x-axis, select <b>Zoom To</b> , type range, then click <b>OK</b> .
Zoom on a specific region of the y-axis	Place the cursor on the left y-axis, then click-drag the Q up or down.
	or
	Right-click the left y-axis, select <b>Zoom To</b> , type maximum, optionally, select <b>Apply to all electropherograms</b> , then click <b>OK</b> .
Unzoom	Double-click the x-axis or y-axis.
	or
	Right-click the x-axis or y-axis, then select <b>Full View</b> .

# Examining Data in the Samples Plot Window

Other tasks you can perform in the Samples Plot window include:

- Adjust the scale of the x-axes (basepairs or data points)
- Adjust the scale of the y-axes (scale to individual maximum, global maximum, or a specific value)
- Show and hide specific dye color peaks
- Display a status line for individual peaks
- Display a Sizing Table, which displays a row of sizing information for each detected peak
- Display a Genotypes Table, which displays a row of genotyping information for each detected peak
- Select peaks, which highlights a corresponding row of data in the Sizing Table

See Figure 2-2 on page 34 for an illustration of some of the above features.

For more information on using the above features, press **F1**, then select the desired topic from the *GeneMapper*<sup>®</sup> *Software Online Help*.

When done viewing the Samples Plot, click X to close the window.

#### **Next Steps**

Create a bin set and generate bins using the Auto Bin feature and reference data as described on page 35.

Chapter 2

Performing the Initial Analysis on the Project

Setting Up the Loss of Heterozygosity Analysis

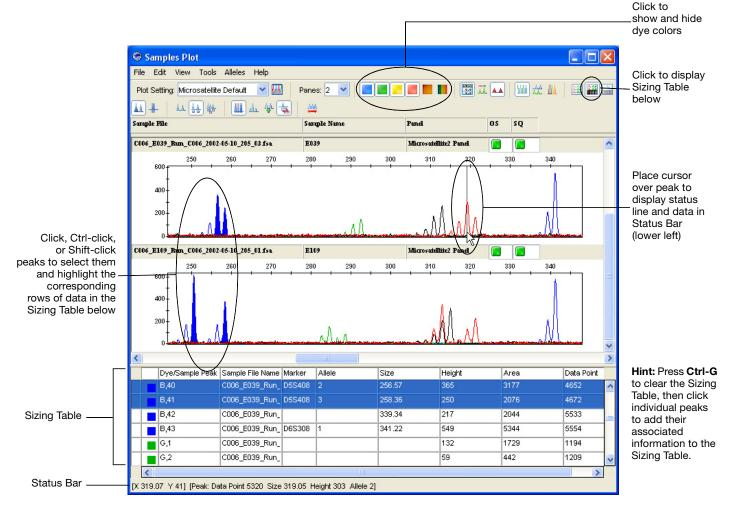


Figure 2-2 Examining and comparing data from different sample files in the Samples Plot

# Creating a Bin Set and Generating Bins Using the Auto Bin Feature

#### Overview

Use the Panel Manager to create bin sets and generate bins (allele definitions).

Before you create a bin set, you must select a kit. You can then associate the bin set with any panels in that kit.

Before you generate bins, you must select a panel and a bin set. Only sample files in projects analyzed with that panel are available to add as reference data to the selected panel to generate the bins. The bins will be associated with markers in the selected panel and stored in the selected bin set.

**Note:** In this guide, you will learn how to create bins using reference data and the Auto Bin feature. However, you can also import bin sets (text files) that contain bin information. Or you can create bins manually. For more information on importing bin sets or creating bins manually, see the *GeneMapper® Software Online Help*.

# Creating a Bin Set

#### To create a bin set:

- 1. Open the Panel Manager by clicking (Tools ▶ Panel Manager).
- **2.** In the Navigation Pane (left), select the **LOH Kit** you created on page 11.
- 3. Click | (Bins ▶ New Bin Set).
- **4.** In the New Bin Set dialog box, type **LOH Bin Set** for the Bin Set Name, then click **OK**.



The LOH Bin Set is added to the Bin Set drop-down list at the top of the Panel Manager. The LOH Bin Set can now be associated with the LOH Panel (or any other panels added to the LOH Kit).

#### Adding Reference Data to a Panel and Bin Set

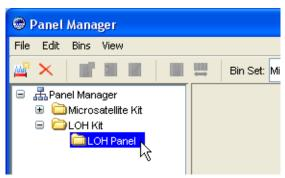
**Note:** You can add all or only a subset of the sample files in a project as reference data. Because the tumor sample files would contain the same alleles as their healthy counterparts, it is not necessary to add all the samples files. In this guide, you will use only the healthy sample files as the reference data for use in creating a panel and bin set.

#### To add reference data to the LOH Panel and LOH Bin Set:

**1.** Make sure the **LOH Bin Set** is selected in the Bin Set drop-down list.



**2.** In the Navigation Pane (left), expand the **LOH Kit**, then select the **LOH Panel** you created on page 11.



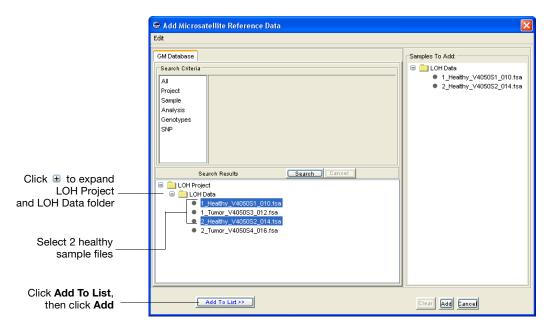
All the markers you created display on the right pane of the Panel Manager (Figure 2-3 on page 38).

#### 3. Click **□** (Bins > Add Reference Data).

The Add Microsatellite Reference Data dialog box opens displaying the LOH Project you created in the lower left pane.

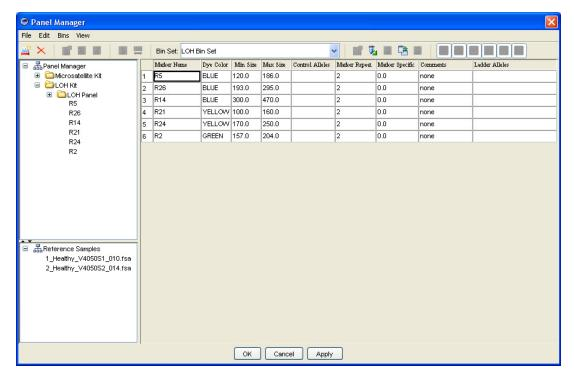
**Note:** The lower left pane always displays all projects that have been analyzed using the selected panel.

- **4.** In the Add Microsatellite Reference Data dialog box:
  - a. Expand LOH Project.
  - **b.** Expand the **LOH** folder.
  - **c.** Press and hold **Ctrl** to select both healthy sample files.
  - d. Click Add to List.
  - e. Click Add.



The selected sample files are added as reference samples to the LOH Panel and appear in the lower half of the Navigation Pane in the Panel Manager (Figure 2-3 on page 38).



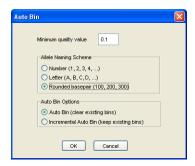


Panel Manager displaying markers and reference Figure 2-3 sample files for selected panel

# Generating Bins using Auto Bin

#### To generate bins using the Auto Bin feature:

- 1. Make sure the LOH Panel is selected in the Navigation Pane and LOH Bin Set is selected in the Bin Set drop-down list, then click (Bins ▶ Auto Bin).
- **2.** In the Auto Bin dialog box, select:
  - Leave Minimum quality value set to **0.1**
  - Rounded basepair for Allele Naming Scheme
  - Auto Bin (clear existing bins) for Auto Bin Options



- 3. Click OK
- **4.** When "Autobinning completed" displays, click **OK**.

#### Reviewing the Markers and Bins

#### To review the markers and bins generated from the reference data:

1. Expand the LOH Panel by clicking 

, then select a marker in the upper half of the Navigation Pane.

A plot (Figure 2-4) displays:

- Marker (pink line)
- Bins (grey columns) for that marker
- Reference alleles (red cross hatches) for each bin
- 2. With a marker selected in the upper half of the Navigation Pane, select a reference sample in the lower half of the Navigation Pane. The plot updates (Figure 2-5) to show the electropherogram peaks for the selected sample.

**Note:** Applied Biosystems recommends selecting each marker to confirm that bins were created for it. If no bins are present, investigate why. See the *GeneMapper® Software Reference and Troubleshooting Guide*.



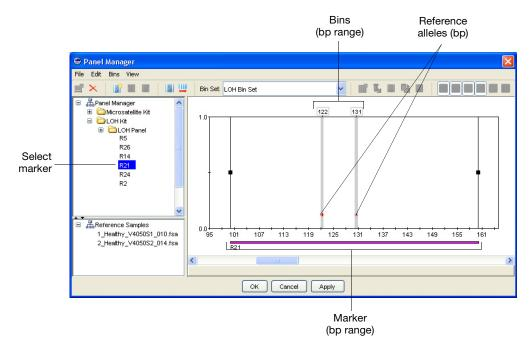


Figure 2-4 Selecting a marker

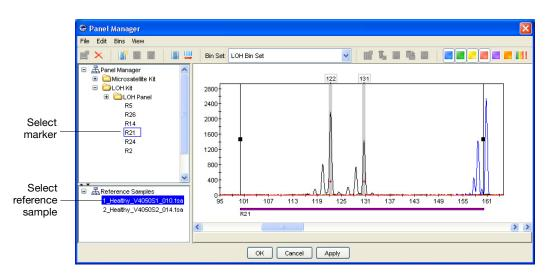


Figure 2-5 Selecting a marker and reference sample

## Accepting the Bin Set

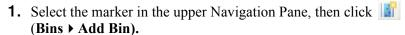
Click **OK** to accept the new bin set and close the Panel Manager.

#### Adding, Editing, and Deleting Bins and Markers (Optional)

To complete the experiment in this guide, you do *not* need to add, edit, or delete any bins or markers. However, you may wish to test these functions by opening the Panel Manager, then selecting the LOH Kit and LOH Panel.

**IMPORTANT!** If you edit or delete any bins or markers, make sure you click **Cancel** at the bottom of the Panel Manager. Clicking OK or Apply can adversely affect the results of the analysis.

#### Adding a Bin to a Marker



- 2. Click in the plot at the location where you want to add the bin.
- **3.** In the Add Bin dialog box, type a **Name**, **Location**, and **Offsets** for the bin, then click **OK**.

#### Editing a Bin

- 1. Click the bin (grey vertical bar) to select it.
- 2. Click (Bins ▶ Edit Bin) or right-click the bin, then select Edit Bin.
- **3.** In the Edit Bin dialog box, edit the **Name**, **Location**, and **Offsets** for the bin, then click **OK**.

#### Editing a Bin Graphically

- **1.** Click the bin (grey vertical bar) to select it (Figure 2-6).
- **2.** Click-drag the blue center line that defines the bin location.
- **3.** Click-drag the left of right handles that define the bin offsets (range).

Note: To correct any undesired change, select Edit ▶ Undo.

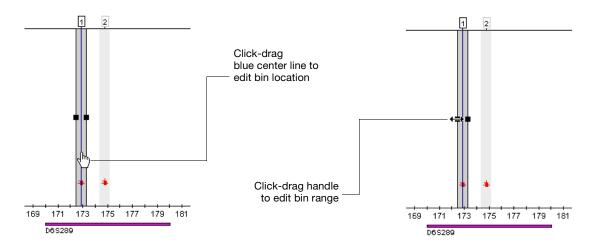


Figure 2-6 Editing a bin graphically

#### Deleting a Bin

To delete a bin from a marker:

- Select the bin, then click (Bins ▶ Delete Bin) or
- Select the bin, right-click the bin, then select **Delete Bin**.

#### Editing a Marker

- **1.** Select the marker in the upper Navigation Pane (Figure 2-7 on page 43).
- **2.** Click-drag the left or right handles that define the marker range.

Note: To correct any undesired change, select Edit ▶ Undo.

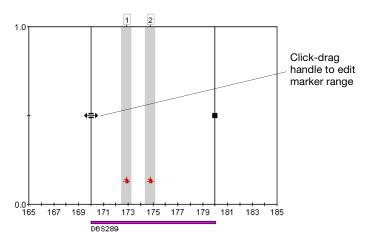


Figure 2-7 Editing a marker

#### Deleting a Marker From a Panel

- **1.** Select the marker in the upper Navigation Pane.
- 2. Click **X** (Edit ▶ Clear Marker).

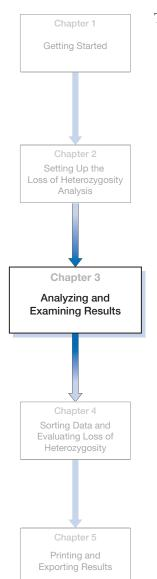
#### **Next Steps**

Analyze and examine the data in the LOH project as described in Chapter 3.

## Chapter 2 Setting Up the Loss of Heterozygosity Analysis Creating a Bin Set and Generating Bins Using the Auto Bin Feature

3

# Analyzing and Examining Results



nis chapter includes:	
Editing the Analysis Method	. 46
Analyzing the Project	. 48
Examining the Results	. 48

## **Editing the Analysis Method**

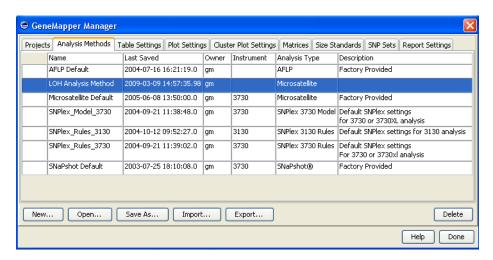
#### Overview

When you created the LOH Analysis Method on page 17, for the initial analysis, you did not select a bin set in the method. Now that you have created a bin set, select that bin set in the analysis method before analyzing the data. The bin set will allow the GeneMapper<sup>®</sup> Software to make allele calls when you analyze the sample files in the project.

#### Editing the Analysis Method

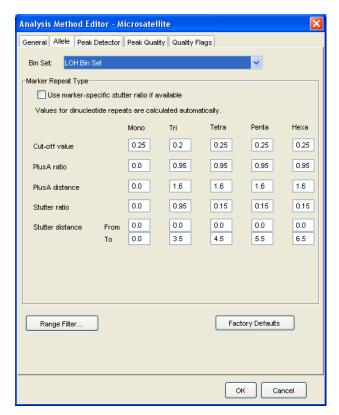
#### To edit the LOH Analysis Method:

- 1. Open the GeneMapper Manager by clicking (Tools ▶ GeneMapper Manager).
- 2. Select the Analysis Methods tab.



3. Select the LOH Analysis Method, then click Open.

**4.** In the Analysis Method Editor, select the **Allele** tab, select **LOH Bin Set** for the Bin Set, then click **OK**.



**5.** Click **Done** to close the GeneMapper Manager.

**Note:** You can also access the Analysis Method Editor by double-clicking any row in the Analysis Method column in the Samples tab of the GeneMapper window (see page 17).

**Next Steps** Analyze the project as described on page 48.

## **Analyzing the Project**

#### Overview

Now that the analysis method specifies a bin set, when you analyze your samples files, the GeneMapper<sup>®</sup> Software will size and genotype the data.

Note the following in the Samples tab of the GeneMapper window:

- The icon displays in the Status column, indicating that the samples are ready to be analyzed and have not been analyzed with the current analysis parameters selected in the Samples tab. This icon displays because you modified the LOH Analysis Method.
- The ✓ icon displays in the REF column, indicating these samples were used as reference data for creating a bin set.

#### Analyzing

Click ► (Analysis ➤ Analyze).

The GeneMapper Software analyzes each sample in the project, displaying its progress in the Status Bar (lower left) of the GeneMapper window.

Next Steps

Examine the results as described below.

## **Examining the Results**

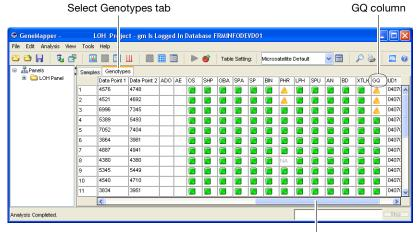
#### Overview

To examine the sizing and genotyping results:

- Review the SQ, associated PQVs, size standard, sample information, and samples plots (described on pages 27 through 33)
- Review the GQ and contributing PQVs (page 49)
- Review the allele calls for each sample (page 50)
- View genotype plots (page 50)
- Examine data in Genotypes Plot window (page 52)
- View project alleles (page 52)

# Reviewing the GQ and PQVs

To review the Genotype Quality (GQ) of the data, select the **Genotypes** tab and scroll to the right.



Click-drag scroll bar to right to view GQ column

If you followed the procedures and used the example data indicated in this guide, the GQ for most samples should be (Pass). The Process Quality Values (PQVs) that contribute the GQ (AN, BD, BIN, LPH, OBA, OS, PHR, SHP, SP, SPA, SPU, and XTLK) should also be , except for a few samples with PHR (Peak Height Ratio). However, yellow PHR is expected for LOH data.

**Hint:** When analyzing LOH data, to reduce the likelihood of PHR being flagged, you may want to edit the Analysis Method to reduce the Minimum Peak Height Ratio in the Heterozygote Balance section of the Peak Quality tab (see page 21).

Investigating Yellow A and Red GQs

**IMPORTANT!** When analyzing your own data, you may find the GQ to be △ (Check) or ♠ (Low Quality) and the contributing PQVs (AN, BD, BIN, LPH, OBA, OS, PHR, SHP, SP, SPA, SPU, and XTLK) to be △, indicating issues with the data, marker or bin definitions, or analysis parameters. To investigate and correct these issues, see the *GeneMapper® Software Reference and Troubleshooting Guide*.

**Note:** Click to sort the samples by GQ score. Samples with a red QQ will be listed at the top of the Genotypes tab.

## Reviewing the Allele Calls

To review the allele calls for each marker in each sample, select the Genotypes tab, then view the Allele 1 and Allele 2 columns.

View allele calls for

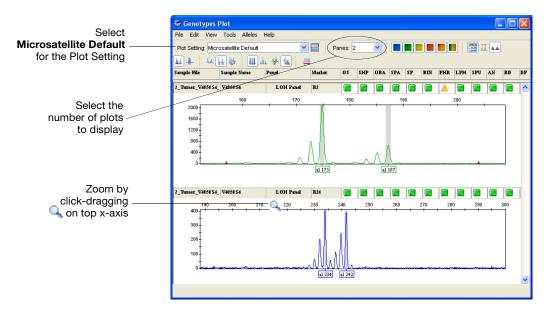
	each mar					marker	•			
Samp	Samples Genotypes									
	Sample File	Sample Name	Panel	Marker	Dye	Allele 1	Allele 2	Size 1	Size 2	Height 1
15	Tumor1_V4050S3_012.fsa	V4050S3	LOH Panel	1Green	G	175	187	174.54	186.95	3014
16	Tumor2_V4050S4_016.fsa	V4050S4	LOH Panel	3Yellow	Y	197	201	196.97	200.96	399
17	Tumor2_V4050S4_016.fsa	V4050S4	LOH Panel	3Blue	В	236	245	236.61	244.89	409
18	Tumor2_V4050S4_016.fsa	V4050S4	LOH Panel	2Yellow	Y	122	131	122.44	130.62	649
19	Tumor2_V4050S4_016.fsa	V4050S4	LOH Panel	2Blue	В	161	161	160.64	160.64	1620
20	Tumor2_V4050S4_016.fsa	V4050S4	LOH Panel	1Yellow	Υ	65	65	64.79	64.79	354
21	Healthy2_V4050S2_014.fsa	V4050S2	LOH Panel	3Blue	В	236	245	236.52	244.86	880
22	Healthy2_V4050S2_014.fsa	V4050S2	LOH Panel	4Blue	В	362	389	361.5	388.71	961
23	Healthy2_V4050S2_014.fsa	V4050S2	LOH Panel	1Green	G	175	187	174.69	187.01	2385
24	Healthy2_V4050S2_014.fsa	V4050S2	LOH Panel	2Yellow	Y	122	131	122.36	130.61	1770

# Viewing Genotype Plots

#### To view the genotypes plots of the samples:

- 1. Select a sample and marker (row) in the Genotypes tab. To select multiple markers, press and hold **Shift** or **Ctrl**. To select all markers, select **Edit** ▶ **Select All**.
- 2. Click (Malysis ➤ Display Plots).

The Genotypes Plot window displays an electropherogram for each selected marker.



#### 3. Select Microsatellite Default for the Plot Setting.

**Note:** Plot Settings control the information displayed in the Genotypes Plot window after analysis. Microsatellite Default is one of the default Plot Settings provided with the GeneMapper Software. You can also edit and create custom Plot Settings in the GeneMapper Manager. For more information, see the *GeneMapper® Software Online Help*.

#### **4.** Zoom on the x- and y-axes in the Genotypes Plot:

То	Then
Zoom on a specific region of the x-axis	Place the cursor on the top x-axis, then click-drag the Q right or left to zoom that plot.
	or
	Right-click the top x-axis, select <b>Zoom To</b> , type range, then click <b>OK</b> .
Zoom on a specific region of the y-axis	Place the cursor on the left y-axis, then click-drag the Q up or down.
	or
	Right-click the left y-axis, select <b>Zoom To</b> , type maximum, optionally, select <b>Apply to all electropherograms</b> , then click <b>OK</b> .
Unzoom	Double-click the x-axis or y-axis.
	or
	Right-click the x-axis or y-axis, then select <b>Full View</b> .

#### Examining Data in the Genotypes Plot Window

Other tasks you can perform in the Genotypes Plot window include:

- Adjust the scale of the x-axes (basepairs or data points)
- Adjust the scale of the y-axes (scale to individual maximum, global maximum, or a specific value)
- Show and hide specific dye color peaks
- Display a status line for individual peaks
- · Add, rename, and delete allele calls
- · Edit and delete markers and bins

For more information on using the above features, press **F1**, then select the desired topic from the *GeneMapper*<sup>®</sup> *Software Online Help*.

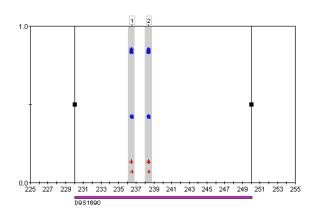
When done viewing the Genotypes Plot, click to close the window.

#### Viewing All Project Alleles

To view all the alleles detected in the sample data for each marker:

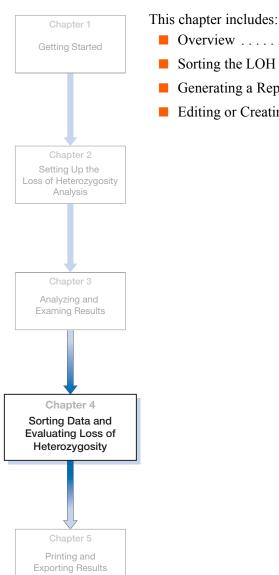
- 1. Open the Panel Manager by clicking ☐ (Tools ▶ Panel Manager).
- **2.** In Navigation Pane, expand the **LOH Kit**, expand the **LOH Panel**, then select a marker in the Navigation Pane (left).
- **3.** Select Bins ▶ Show Project Alleles.

The project alleles (alleles detected in the sample data) appear as blue asterisks \* in each bin. The y-axis position of each \* indicates the GQ score for that marker and sample.





# Sorting Data and Evaluating Loss of Heterozygosity



1	
Overview	54
Sorting the LOH Data	56
Generating a Report to Calculate and Evaluate LOH	58
Editing or Creating a Custom LOH Report Setting	59

### **Overview**

An LOH analysis investigates and compares healthy tissue and suspected diseased tissue from the same individual. Both the healthy and diseased samples should show two alleles for a given heterozygous microsatellite marker. If LOH is occurring in the diseased tissue, it will be evident by a decrease in peak height of one of the two alleles (relative to the healthy allele peak heights). The reason the tumor sample shows a decrease in peak height (instead of an absence of the peak) is that during isolation, the tumor sample is contaminated with healthy cells, thus introducing some wild-type DNA into the tumor specimen.

The Report Manager in the GeneMapper Software includes an LOH Default report setting that calculates and compares peak height ratios of microsatellite alleles between healthy and diseased tissues (Figure 4-1). It then evaluates and flags any samples that show potential LOH based on set peak-height ratio thresholds, so these samples can be further investigated. You can alter these thresholds based on the observed level of wild-type DNA contamination in the tumor samples.

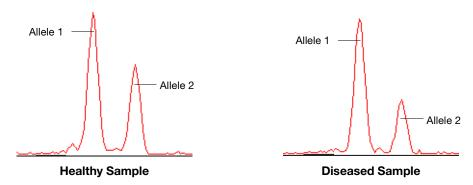


Figure 4-1 Diseased sample shows loss in peak height of Allele 2

The calculations and LOH criteria used by the LOH Default report setting are summarized below:

 Calculates the Allele Ratio between 2 allele peaks for each marker for each sample using the following equation:

Allele Ratio = 
$$\frac{\text{Peak Height of Allele 1}}{\text{Peak Height of Allele 2}}$$

Calculates Allelic Imbalance between Allele Ratios for the healthy sample and diseased sample for each marker using the following equation:

3. Evaluates the Allelic Imbalance for each marker and sample pair then flags those with scores above or below set thresholds:

If Allelic Imbalance > 1.35, then flag as LOH Candidate
OR
If Allelic Imbalance < 0.67, then flag as LOH Candidate

Figure 4-2 Calculations and thresholds used by the LOH Default report setting

**IMPORTANT!** You can edit the LOH Default report setting or create your own report setting as appropriate for your data (see page 59).

## Sorting the LOH Data

#### Overview

Before you can apply the LOH Default report setting to calculate and evaluate potential LOH, it is critical to sort the data so that healthy samples and diseased samples for each marker are listed consecutively.

#### Sorting To sort the LOH data:

- 1. Select the **Genotypes tab** in the GeneMapper window.
- **2.** Sort by sample files by pressing and holding the **Shift** key, then clicking the Sample File column header.

All markers from the 1 Healthy sample file are listed first, followed by all samples from the 1 Tumor sample file, and so on.

		Shift-click
	Sample File ——	Sample File
1	1_Healthy_V4050S1_010.fsa	header
2	1_Healthy_V4050S1_010.fsa	
3	1_Healthy_V4050S1_010.fsa	
4	1_Healthy_V4050S1_010.fsa	
5	1_Healthy_V4050S1_010.fsa	
6	1_Healthy_V4050S1_010.fsa	
7	1_Tumor_V4050S3_012.fsa	
8	1_Tumor_V4050S3_012.fsa	
9	1_Tumor_V4050S3_012.fsa	
10	1_Tumor_V4050S3_012.fsa	
11	1_Tumor_V4050S3_012.fsa	
12	1_Tumor_V4050S3_012.fsa	
13	2_Healthy_V4050S2_014.fsa	
14	2_Healthy_V4050S2_014.fsa	
15	2_Healthy_V4050S2_014.fsa	
16	2_Healthy_V4050S2_014.fsa	

Shift aliak

Figure 4-3 Sorting by sample file

**3.** Sort by Marker by pressing and holding the **Shift** key, then clicking the Marker column header.

This sorts the results so that healthy samples and diseased samples for the same marker are listed consecutively.

					Shift-click
	Sample File	Sample Name	Panel	Marker —	Marker
1	1_Healthy_V4050S1_010.fsa	V4050S1	M2 LOH Panel	R14	header
2	1_Tumor_V4050S3_012.fsa	V4050S3	M2 LOH Panel	R14	
3	2_Healthy_V4050S2_014.fsa	V4050S2	M2 LOH Panel	R14	
4	2_Tumor_V4050S4_016.fsa	V4050S4	M2 LOH Panel	R14	
5	1_Healthy_V4050S1_010.fsa	V4050S1	M2 LOH Panel	R2	
6	1_Tumor_V4050S3_012.fsa	V4050S3	M2 LOH Panel	R2	
7	2_Healthy_V4050S2_014.fsa	V4050S2	M2 LOH Panel	R2	
8	2_Tumor_V4050S4_016.fsa	V4050S4	M2 LOH Panel	R2	
9	1_Healthy_V4050S1_010.fsa	V4050S1	M2 LOH Panel	R21	
10	1_Tumor_V4050S3_012.fsa	V4050S3	M2 LOH Panel	R21	
11	2_Healthy_V4050S2_014.fsa	V4050S2	M2 LOH Panel	R21	
12	2_Tumor_V4050S4_016.fsa	V4050S4	M2 LOH Panel	R21	
13	1_Healthy_V4050S1_010.fsa	V4050S1	M2 LOH Panel	R24	
14	1_Tumor_V4050S3_012.fsa	V4050S3	M2 LOH Panel	R24	1
15	2_Healthy_V4050S2_014.fsa	V4050S2	M2 LOH Panel	R24	1
16	2_Tumor_V4050S4_016.fsa	V4050S4	M2 LOH Panel	R24	

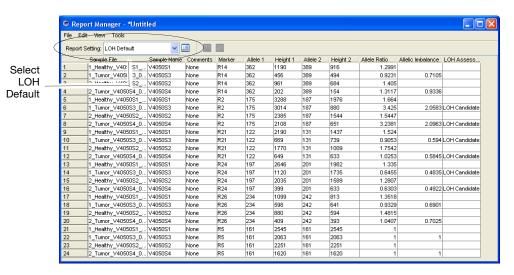
Figure 4-4 Sorting by sample file and marker

**Next Steps** Generate a report to calculate and evaluate the LOH as described on page 58.

## Generating a Report to Calculate and Evaluate LOH

To generate a report that calculates and evaluates loss of heterozygosity:

- **1.** Make sure you sorted the data as described on page 56.
- 2. Select all the samples in the Genotypes tab by selecting Edit ➤ Select All.
- 3. Select Analysis ➤ Report Manager.
- **4.** In the Report Manager dialog box, select **LOH Default** from the Report Setting drop-down list.



The LOH Default report setting is applied to the LOH Project data. The resulting report shows the Allele Ratio, Allelic Imbalance, and LOH Assessment for each sample or pair of samples.

**5.** Save, print, or export the generated report by selecting the appropriate command from the File menu.

**IMPORTANT!** When generating a report for your own data, make sure samples do not have a Low Quality SQ score in the Samples tab. If they do, the samples and associated markers appear in the Genotypes tab, but with no data to be used by the report setting.

## **Editing or Creating a Custom LOH Report Setting**

#### Overview

The LOH Default report setting provided with the GeneMapper Software assumes the following:

- Allelic Imbalance Row Calculation When calculating the Allelic Imbalance (ratio of peak-height ratios) between two samples, the report setting compares every two rows in the Genotypes tab. This calculation assumes your sample files are named so that healthy samples and diseased samples for each individual and marker are listed consecutively (in pairs) after being sorted by sample file and marker.
- **LOH Assessment Thresholds** The threshold values used to flag samples as LOH candidates are > 1.35 and < 0.67.

You can edit the above settings in the LOH Default report setting as appropriate for your samples and experiment.

#### Editing the Allelic Imbalance Row Calculation

Depending on how your sample files are named, healthy and tumor files from the same individual may or may not list consecutively after being sorted by sample file and marker (Figure 4-5).

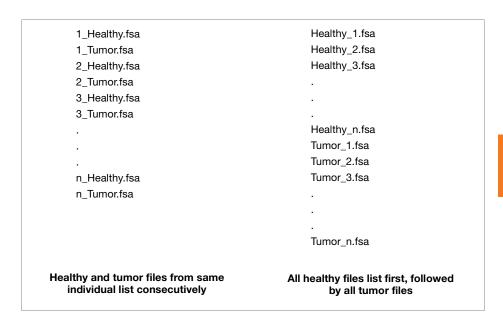


Figure 4-5 How sample file naming convention affects the order files list after being sorted by sample and marker

#### To change the row calculation for the Allelic Imbalance in the LOH Default report setting:

- **1.** Open the GeneMapper Manager by clicking (Tools ▶ GeneMapper Manager).
- **2.** Select the **Report Settings** tab (Figure 4-5).
- 3. Select the LOH Default report setting, then click Open.

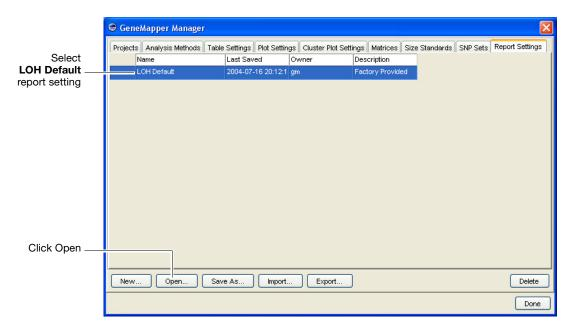


Figure 4-6 GeneMapper Manager Report Settings tab

- **4.** In the Report Settings Editor (Figure 4-5):
  - a. Select the Calculation tab.
  - b. Select Allelic Imbalance from the Selected Columns list (right pane).
  - **c.** Click to move Allelic Imbalance to the Calculation tab.

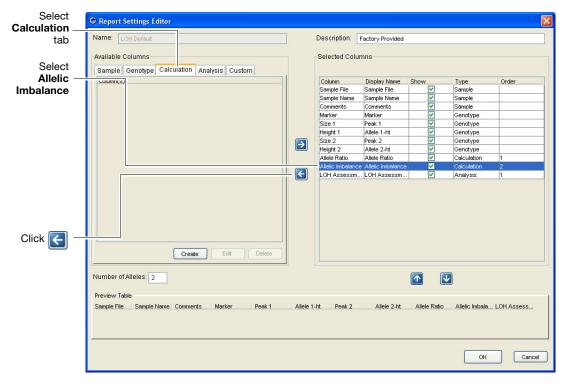


Figure 4-7 Report Settings Editor

- 5. Select Allelic Imbalance in the Calculation tab, then click Edit.
- **6.** In the Edit Calculation dialog box, replace 2 with n + 1 in the two fields shown in Figure 4-5.

**Note:** The value *n* is the sample set number, that is the number of healthy samples or the number of tumor samples (see Figure 4-5 on page 59). For example, if your sample set contains 5 healthy and 5 tumor samples, you would type 6.



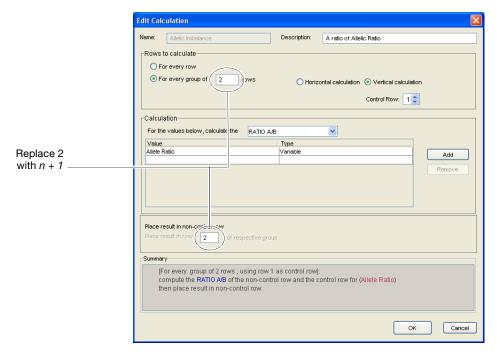


Figure 4-8 Edit Calculation dialog box

- 7. Click **OK** to close the Edit Calculation dialog box.
- **8.** In the Report Settings Editor (Figure 4-5), click to move Allelic Imbalance from the Calculation tab to the Selected Columns list.
- **9.** In the Report Settings Editor (Figure 4-5):
  - a. Select the Analysis tab.
  - **b.** Select **LOH Assessment** from the Selected Columns list (right pane).
  - **c.** Click to move LOH Assessment to the Analysis tab.
- 10. Select LOH Assessment in the Analysis tab, then click Edit.

**11.** In the Analysis section of the Edit Analysis dialog box, select **For every group of**, then type n + 1 in the two fields shown in Figure 4-5.

**Note:** The value *n* is the sample set number, that is the number of healthy samples or the number of tumor samples (see Figure 4-5 on page 59). For example, if your sample set contains 5 healthy and 5 tumor samples, you would type 6.

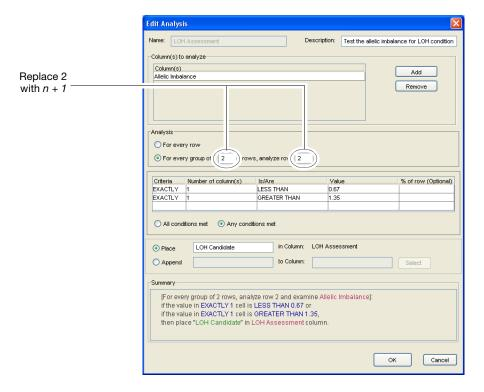


Figure 4-9 Edit Analysis dialog box

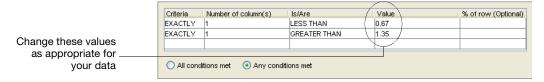
- **12.** Click **OK** to close the Edit Analysis dialog box.
- **13.** In the Report Settings Editor (Figure 4-5 on page 61), click to move LOH Assessment from the Analysis tab to the Selected Columns list.
- **14.** Click **OK** to close the Report Settings Editor. Click **Done** to close the GeneMapper Manager.

#### Editing the LOH Assessment Thresholds

Select LOH Assessment thresholds appropriate for the observed level of wild-type DNA contamination in tumor samples. For more information, see the discussion on page 54.

To change the thresholds for the LOH Assessment in the LOH Default report setting:

- 1. Open the GeneMapper Manager by clicking (Tools ▶ GeneMapper Manager).
- **2.** Select the **Report Settings** tab (Figure 4-5 on page 60).
- **3.** Select the **LOH Default** report setting, then click **Open**.
- **4.** In the Report Settings Editor (Figure 4-5 on page 61):
  - a. Select the Analysis tab.
  - **b.** Select **LOH Assessment** from the Selected Columns list (right pane).
  - **c.** Click to move LOH Assessment to the Analysis tab.
- **5.** Select LOH Assessment in the Analysis tab, then click **Edit**.
- **6.** In the Edit Analysis dialog box, in the Value column, replace the lower and upper thresholds of 0.67 and 1.35 with thresholds appropriate for your data.



- 7. Click **OK** to close the Edit Analysis dialog box.
- **8.** In the Report Settings Editor (Figure 4-5), click to move LOH Assessment from the Analysis tab to the Selected Columns list.
- **9.** Click **OK** to close the Report Settings Editor. Click **Done** to close the GeneMapper Manager.

5

## **Printing and Exporting Results**

Chapter 1			
Getting Started			
Chapter 2			
Setting Up the			
Loss of Heterozygosity  Analysis			
Aildiyələ			
Chapter 3			
Analyzing and			
Examing Results			
Chapter 4			
Sorting Data and			
Evaluating Loss of			
Heterozygosity			
Chapter 5			
Printing and			
Exporting Results			

This chapter	includes:
--------------	-----------

### **Printing Results**

You can print results from the following windows and tabs by selecting **File > Print**:

Window/Tab	Access From GeneMapper Project Window by selecting
GeneMapper window – Samples tab	View ▶ Samples
GeneMapper window – Genotypes tab	View ▶ Genotypes
GeneMapper window – Info tab	View ▶ Sample Info
GeneMapper window – Raw Data tab	View ▶ Raw Data
GeneMapper window – EPT Data tab	View ▶ EPT Data
Samples Plot window	The Samples tab, then <b>Analysis ▶ Display</b> Plots
Genotypes Plot window	The Genotypes tab, then Analysis ➤ Display Plots
Report Manager window	Analysis ▶ Report Manager

**Note:** You can also print reports. For information on creating report settings and generating reports, see the  $GeneMapper^{\otimes}$  Software  $Online\ Help$ .

### **Exporting Results**

#### Exporting Samples Tab and Genotypes Tab

To export the results displayed in the Samples tab and Genotypes tab of the GeneMapper window:

- **1.** Prepare the content and format of the data to export:
  - a. Select the desired Table Setting from the drop-down list at the top of the GeneMapper window. The Table Setting controls which columns display and the sorting order for the samples.
  - b. Optionally, sort the data to determine the order that the samples appear. Select Edit ➤ Sort or Shift-click the column header in the Samples tab or Genotypes tab. You can also click (Analysis ➤ Low Quality on Top) to sort the samples by GQ score.

**Note:** For more information on editing or creating Table Settings and sorting data, see the *GeneMapper*<sup>®</sup> *Software Online Help*.

- **2.** Select one of the following commands:
  - File > Export Table Exports information displayed in the selected tab.
  - File > Export Combined Table Exports information displayed in both tabs. (This command is available only when the Samples tab is selected.)

#### Exporting Kits To

To export all kits in the Panel Manager:

- 1. Open the Panel Manager by clicking ☐ (Tools ▶ Panel Manager).
- 2. Select File > Export All Kits.

#### **Exporting Panels**

To export all panels in a kit:

- 1. Open the Panel Manager by clicking (Tools ▶ Panel Manager).
- **2.** Select the kit in the Navigation Pane (left).
- **3.** Select File ▶ Export Panels.

## Exporting Bin Sets

#### To export a bin set:

- 1. Open the Panel Manager by clicking ☐ (Tools ▶ Panel Manager).
- **2.** In the Navigation Pane, select the kit with which the bin set is associated.
- **3.** Select the bin set from the Bin Set drop-down list.
- 4. Select File ▶ Export Bin Set.

#### Exporting Projects, Methods, Settings, and Size Standards

To export projects, analysis methods, table settings, plot settings, reports settings, and size standards:

- 1. Open the GeneMapper Manager by clicking ☐ (Tools ▶ GeneMapper Manager).
- **2.** Select one of the following tabs:
  - Projects
  - · Analysis Methods
  - · Table Settings
  - Plot Settings
  - Report Settings
  - Size Standards
- **3.** Select the object(s) you want to export. Press and hold **Shift** or **Ctrl** to select multiple objects.
- 4. Click Export.

# Exporting Reports

You can also export reports. For information on creating report settings and generating reports, see the *GeneMapper® Software Online Help*.

### Index

A	Auto Bin 39
adding bins to markers 41 sample files to project 15 size standard labels 30	B bin adding to a marker 41
Advanced peak detection algorithm 20 allele calls, reviewing 50 Allele Ratio 55 Allele tab of Analysis Method Editor 19, 47 alleles, project 52 Allelic Imbalance 55, 59 analysis method creating 17 editing 46 exporting 68 saving 23 selecting bin set 19, 47	creating manually 35 definition 7, 35 deleting 42 editing 41, 42 generating using Auto Bin 39 reviewing 39 bin set accepting 41 creating 35 definition 7, 35 exporting 68 importing 35 selecting in analysis method 19, 47
Analysis Method Editor Allele tab 19, 47	bold text, when to use v
General tab 18 Peak Detector tab 20 Peak Quality tab 21 Quality Flags tab 22 analysis parameters definition 7, 17 setting 17 analysis. See LOH microsatellite analysis	conventions bold text v for describing menu commands v IMPORTANTS! vi in this guide v italic text v Notes vi
analyzing a project 26, 48	user attention words vi
Applied Biosystems contacting viii customer feedback on documentation viii Information Development department viii Technical Support viii assumptions for using this guide v	creating analysis method 17 bin set 35 bins 39 kit 11 markers 11 panel 11 project 15

report setting 68 customer feedback, on Applied Biosystems documents viii	F files. See sample files fill down 24
D	G
data	G
See example data	General tab of Analysis Method Editor 18
See reference data	generating
See sample files	report for LOH 58
deleting	reports 68
bins 42	genotype quality. See GQ
markers 43	Genotypes Plot window
size standard labels 30	displaying 50
disclaimer, license ii	examining data in 52
documentation, related vii	exporting 67
documentation, related vii	printing 66
_	zooming 51
E	Genotypes tab 49
editing	Genotypes Table 33
Allelic Imbalance 59	GQ
analysis method 46	investigating 49
bins 41, 42	reviewing 49
LOH Assessment Thresholds 64	Teviewing 19
markers 43	Н
report setting 59	П
size standard labels 30	help, online, accessing vii, viii
example data	hiding dye color peaks 33, 52
instrument used 4	
marker information 5, 14	1
overview 4	•
size standard for 4	importing
exporting	bin set 35
analysis methods 68	panels 11
bin sets 68	Information Development department,
Genotypes Plot window 67 kits 67	contacting viii
panels 67	instruments
plot settings 68	compatible with LOH microsatellite
projects 68	analyses 3
report settings 68	for example data 4
reports 58, 68	italic text, when to use v
results 67	,
Samples Plot window 67	K
size standards 68	
table settings 68	kits

creating 11 definition 7, 11	overriding SQ 30
exporting 67 type 12	Р
L	Panel Manager opening 11
license disclaimer ii	panels
	adding reference data 36
LOH definition 2, 54 evaluating 53	creating 11 definition 7, 11 exporting 67
LOH Assessment thresholds 55, 59, 64	importing 11
LOH Default report setting 55, 59	selecting for analysis 23
LOH microsatellite analysis	peak detection algorithm 20
compatible instruments 3 definition 2, 54	Peak Detector tab of Analysis Method Editor 20
evaluating LOH 53 flowchart 6	Peak Quality tab of Analysis Method Editor 21
sorting data 56	plot setting 31, 51, 68
	Plot window
M markers	See Genotypes Plot window See Samples Plot window
adding bins to 41	plots. See sample plots
creating 11	primers, custom 3
definition 2, 7, 11	printing
deleting from a panel 43 editing 43 for example data 5, 14 plots, viewing 50 reviewing 39 viewing plots 50	Genotypes Plot window 66 reports 58, 66 results 66 Samples Plot window 66 project
menu commands, conventions for describing v	adding sample files to 15 analyzing 26, 48
method. See analysis method	creating 15
-	exporting 68
microsatellite analysis definition 2 markers 2	saving 25 setting analysis parameters 17 setting table settings 17
See also LOH microsatellite analysis setting up 9	project alleles viewing 52
MSDSs, obtaining viii	5
	Q
0	Quality Flags tab of Analysis Method
online help, accessing vii, viii	Editor 22

R	examining data in 33
reference data	exporting 67
adding to a panel 36	printing 66 zooming 32
icon 48	Samples tab 17
Report Manager 58	saving
report setting	analysis method 23
creating 68	project 25
editing 59	reports 58
LOH Default 55, 59	setting up analysis 9
Report Settings Editor 60	showing and hiding dye color peaks 33, 52
report settings, exporting 68	size calling curve 29
report variables	Size Match Editor 28
Allele Ratio 55	size quality. See SQ
Allelic Imbalance 55, 59	
LOH Assessment 55, 59, 64	size standard
reports	custom 30 examining 28
exporting 58, 68 generating 68	exporting 68
generating for LOH 58	for example data 4
printing 58, 66	labels, adding 30
saving 58	labels, deleting 30
results	labels, editing 30
exporting 67	selecting for analysis 23 troubleshooting 30
printing 66	Sizing Table 33
S	software starting and logging in 7
sample files	terms defined 7
adding to project 15	sorting
location 4	LOH samples 56
naming convention 4, 59	samples 27, 49
plots, viewing 31, 50	SQ
sorting 27, 49, 56 viewing information on 30	investigating 27
viewing plots 31, 50	overriding 30
zooming 32, 51	reviewing 27
sample plots	score 28
viewing 31, 50	Status Bar 26
zooming 32, 51	
samples	Т
sorting 27, 49	table settings 17, 24, 68
sorting for LOH evaluation 56	Technical Support, contacting viii
zooming 32, 51	text conventions v
Samples Plot window	thresholds. See LOH Assessment thresholds
	unesholds, see LOH Assessment unesholds

```
training, information on viii
troubleshooting the size standard 30
U
unzooming 32, 51
user attention words, described vi
X
x-axis
   scale 33, 52
   zooming 32, 51
Y
y-axis
   scale 33, 52
   zooming 32, 51
Z
zooming
   Genotypes Plot window 51
   Samples Plot window 32
   unzooming 32, 51
   x-axis 32, 51
   y-axis 32, 51
```

Index

