

Microsatellite Analysis Getting Started Guide





Getting Started

2

Setting Up the Microsatellite Analysis

3

Analyzing and Examining Results



Printing and Exporting Results

2



Microsatellite Analysis Getting Started Guide

Getting Started

Setting Up the Microsatellite Analysis

Analyzing and Examining

Results

3

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

Information in this document is subject to change without notice.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. TO THE FULLEST EXTENT ALLOWED BY LAW, IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF, WHETHER OR NOT FORESEEABLE AND WHETHER OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

NOTICE TO PURCHASER: DISCLAIMER OF LICENSE

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.

The AFLP process is covered by patents owned by Keygene N.V.

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.

AFLP is a registered trademark of Keygene N.V.

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.

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 4403672 Rev. A 04/2009

Contents

	Preface	V
	How to Use This Guide	٠
	How to Obtain More Information	v
	How to Obtain Support	vii
Chapter 1	Getting Started	1
	About Microsatellite Analyses	2
	About the Example Data	
	Microsatellite Analysis Workflow	4
	GeneMapper® Software Terms	5
	Starting the Software and Logging In	5
	Using This Guide With Your Own Sample Files	6
	Alternatives to the Procedures in This Guide	7
Chapter 2	Setting Up the Microsatellite Analysis	g
	Overview	10
	Creating a Kit, Panel, and Markers	
	Creating a New Project and Adding Sample Files	
	3 , 3 ,	
	Setting Analysis Parameters and Table Settings for the Project	18
	Setting Analysis Parameters and Table Settings for the Project Performing the Initial Analysis on the Project	
	Performing the Initial Analysis on the Project	27
	Performing the Initial Analysis on the Project	27
Chapter 3	Performing the Initial Analysis on the Project	27
Chapter 3	Performing the Initial Analysis on the Project	36 36

Contents

	Analyzing the Project	
Chapter 4	Printing and Exporting Results	53
	Printing Results	54
	Exporting Results	55
	Index	57

Preface

How to Use This Guide

Purpose of This Guide

The GeneMapper® Software Version 4.1 Microsatellite Analysis Getting Started Guide provides brief, step-by-step instructions for sizing and genotyping 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® XP 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 DVD*.

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 http://www.appliedbiosystems.com, then click the link for Support.

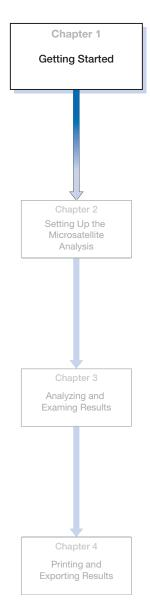
At the Support page, you can:

- 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



About Microsatellite Analyses	2
About the Example Data	2
■ Microsatellite Analysis Workflow	4
■ GeneMapper® Software Terms	5
■ Starting the Software and Logging In	5
■ Using This Guide With Your Own Sample Files	6
■ Alternatives to the Procedures in This Guide	7

This chapter includes:

About 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.

Custom Primers

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

Compatible Kits and Instruments

For information about Applied Biosystems chemistry kits and electrophoresis instruments that are compatible with microsatellite analyses, see the *GeneMapper® Software Version 4.1 Quick Reference Guide* (PN 4403615).

About the Example Data

Sample File Location

To perform the exercise described in this guide, use the three sample files (.fsa) located on your computer at:

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

Note: The above location will vary depending on the installation of the GeneMapper[®] Software. The default installation is the D drive.

Instrument and Size Standard

Sample files were generated by running PCR-amplified and fluorescently tagged microsatellite samples on an ABI PRISM[®] 3100 Genetic Analyzer using the GeneScan[™] 500 LIZ[®] Size Standard.

Marker Information

The sample files contain the following 17 fluorescently tagged markers.

Marker	Dye Label	Allele Size Range (bp)
D6S264	Blue	108 – 130
D6S1574	Blue	146 – 172
D6S276	Blue	201 – 233
D5S408	Blue	240 – 285
D6S308	Blue	326 – 354
D6S287	Green	104 – 138
D6S292	Green	154 – 176
D6S434	Green	201 – 245
D5S426	Green	274 – 298
D5S1981	Yellow	112 – 125
D6S257	Yellow	167 – 195
D6S446	Yellow	212 – 234
D5S641	Yellow	299 – 339
D5S433	Red	69 – 99
D5S406	Red	168 – 196
D5S400	Red	221 – 243
D6S309	Red	307 – 333

Note: You will use the information in the above table when creating a panel and markers in Chapter 2, "Setting Up the Microsatellite Analysis."

Microsatellite Analysis Workflow

The following flowchart summarizes the steps for performing a microsatellite analysis using the GeneMapper® Software:

Set Up the Microsatellite 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.
- 5. 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.

Print and Export the Results (Optional) (Chapter 4)

- · 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 [®] 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 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 data using the GeneMapper[®] 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

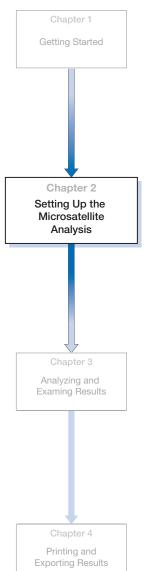
The GeneMapper Software includes an Autoanalysis feature that can eliminate most of the tasks leading up to the analysis of a microsatellite project. Much of Chapter 2, "Setting Up the Microsatellite Analysis," explains how to manually create, add samples to, and analyze projects for use in microsatellite 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 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 projects because it automate many of the tasks explained in Chapter 2, "Setting Up the Microsatellite 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 Microsatellite 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 f or the Project	18
	Performing the Initial Analysis on the Project	27
	Creating a Bin Set and Generating Bins Using the Auto Bin Feature	36

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 \(\overline{\overli pressing F1.



Creating a Kit, Panel, and Markers

Overview

You create the following hierarchical objects in the Panel Manager:

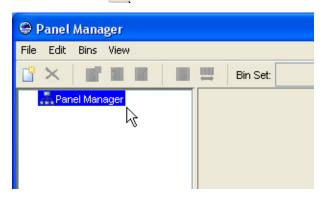
- **Kit** A group of panels
- Panel A group of markers
- Marker A fragment size range (bp), dye color, and repeat length

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[®] Software for some of the LMS kits available from Applied Biosystems. For more information on importing panels, see the *GeneMapper*[®] *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).

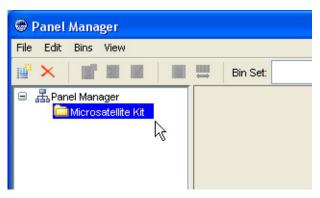


3. In the New Kit dialog box, type **Microsatellite Kit** for the Kit Name, select **Microsatellite** for the Kit Type, then click **OK**.



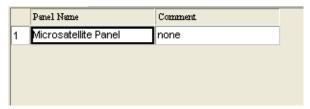


Microsatellite Kit appears in the Navigation Pane (left side).



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

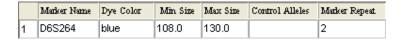
5. In the right pane of the Panel Manager, select **New Panel**, type **Microsatellite Panel** for the Panel Name, then press **Enter**.



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



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



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

Marker	Dye Color	Minimum Size (bp)	Maximum Size (bp)	Marker Repeat
D6S1574	Blue	146	172	2
D6S276	Blue	201	233	2
D5S408	Blue	240	285	2
D6S308	Blue	326	354	2
D6S287	Green	104	138	2
D6S292	Green	154	176	2
D6S434	Green	201	245	2
D5S426	Green	274	298	2
D5S1981	Yellow	112	125	2
D6S257	Yellow	167	195	2
D6S446	Yellow	212	234	2
D5S641	Yellow	299	339	2
D5S433	Red	69	99	2
D5S406	Red	168	196	2
D5S400	Red	221	243	2
D6S309	Red	307	333	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.

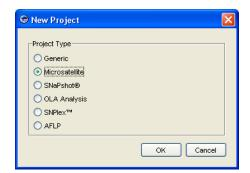
Creating a New Project and Adding Sample Files

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).



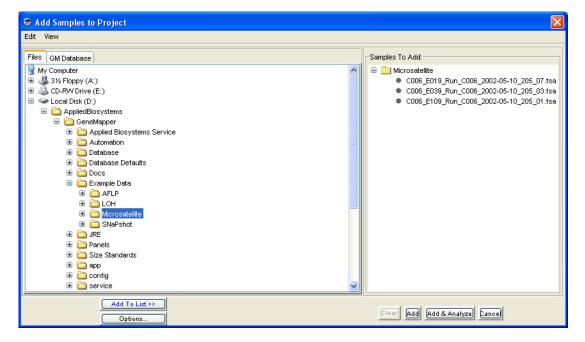
- 2. In the New Project dialog box, select **Microsatellite**, then click **OK**
- **4.** In the Add Samples to Project dialog box, in the Files tab, navigate to:

 $\label{lem:decomp} $$ \down{10pt} $$$ \down{10pt} $$$$ \down{10pt} $$$ \down{10pt} $$$$ \$

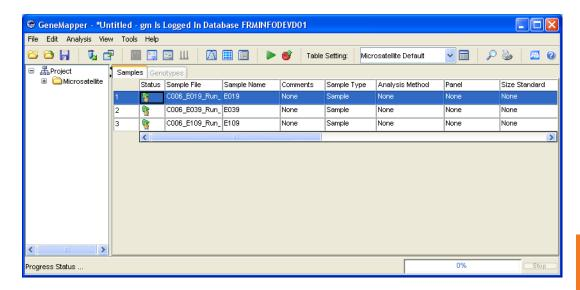
Note: The above location will vary depending on the installation of the GeneMapper[®] Software. The default installation is the D drive.

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

Note: For this guide you added all three sample files in the Microsatellite 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 three sample files from the Microsatellite 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 18.

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

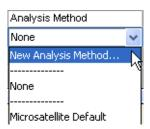
You set analysis parameters that determine the peak detection, sizing, and genotyping algorithms used by the GeneMapper[®] Software to analyze all sample files in a project.

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.

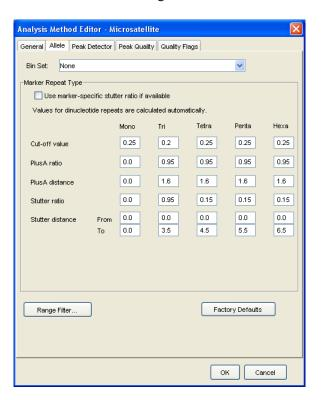
3. In the New Analysis Method dialog box, select **Microsatellite** for Analysis Type, then click **OK**.



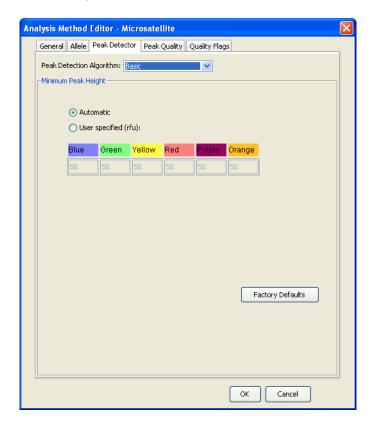
- **4.** In the Analysis Method Editor dialog box, select and edit the five tabs.
 - **General** This tab includes reference information about the method. Type **Microsatellite Analysis Method** for the Name. Optionally, type a description and the instrument on which the data was generated.



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

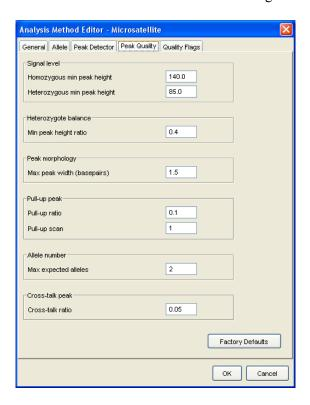


• **Peak Detector** – This tab includes settings that determine peak detection and sizing of peaks. Select **Basic** for the Peak Detection Algorithm. Leave the default values for all other settings.



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

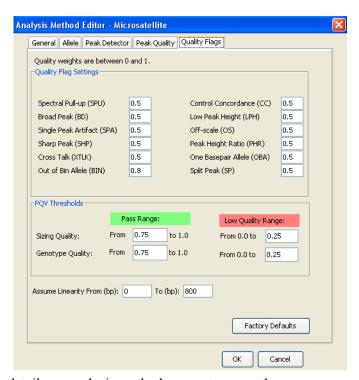
Type **140.0** for the Homozygous min peak height. Type **85.0** for the Heterozygous min peak height. Type **0.4**. for the Min peak height ratio. Leave the default values for all other settings.



• 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).

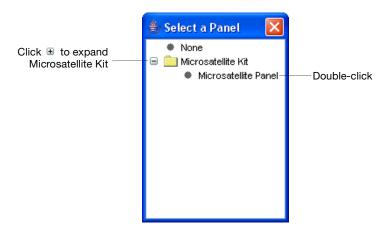
Leave the default values for all settings.



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



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



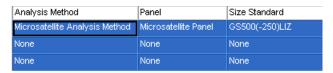
3. Select the first row in the Size Standard column, then select GS500(-250)LIZ from the drop-down list.

Note: In the GeneMapper software, the following size standards are available for use with samples run with the GeneScan[™] 500 LIZ® Size Standard:

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

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.

- **4.** 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.



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.

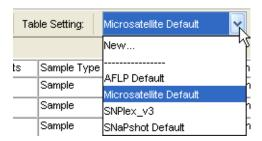


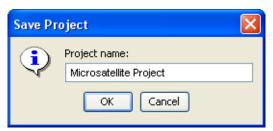
Table Settings control the information displayed in the Samples tab and Genotypes tab after analysis. Microsatellite Default is one of the default Table Settings provided with the GeneMapper Software.

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 **Microsatellite Project**, then click **OK**.



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

Next Steps

Perform the initial analysis on the project as described on page 27.

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[®] 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

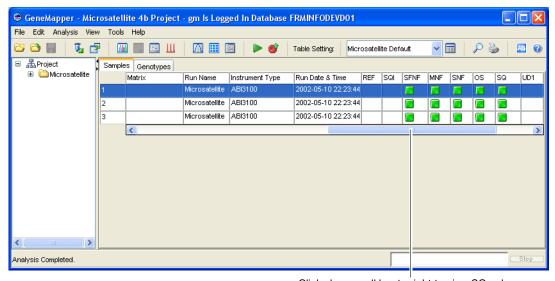
Click ► (Analysis ➤ Analyze).

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). The Process Quality Values (PQVs) that contribute to the SQ (SFNF, MNF, SNF, and OS) should also be .

Investigating Yellow 🛕 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 29.

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.



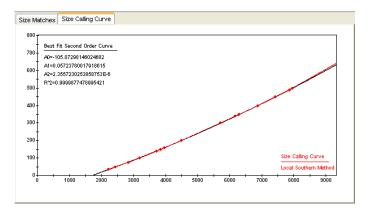
Figure 2-1 Size Match Editor - Size Matches tab

- **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 31.

- **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 31.
- **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.

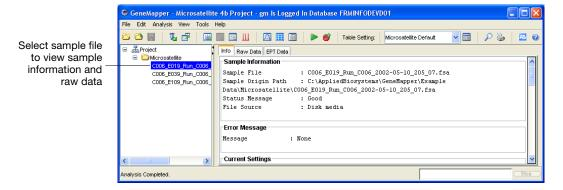
Table 2-1 Troubleshooting the size standard

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 Override SQ 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 Apply to reanalyze the data with the updated sizing information. For more information, see the <i>GeneMapper</i> ® <i>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 $GeneMapper^{\mathbb{R}}$ 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

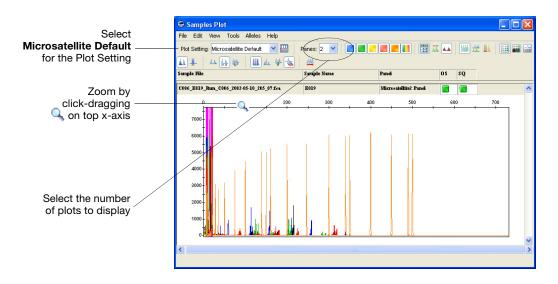


Viewing Sample Plots

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 Shift while click-dragging to zoom only the selected plot.
	or
	Right-click the top x-axis, select Zoom To , type range, then click OK .
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 Zoom To , type maximum, optionally, select Apply to all electropherograms , then click OK .
Unzoom	Double-click the x-axis or y-axis.
	or
	Right-click the x-axis or y-axis, then select Full View .

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 35 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*[®] *Software Online Help*.

When done viewing the Samples Plot, click \(\subseteq \text{to close the window.} \)

Next Steps

Create a bin set and generate bins using the Auto Bin feature and reference data as described on page 36.

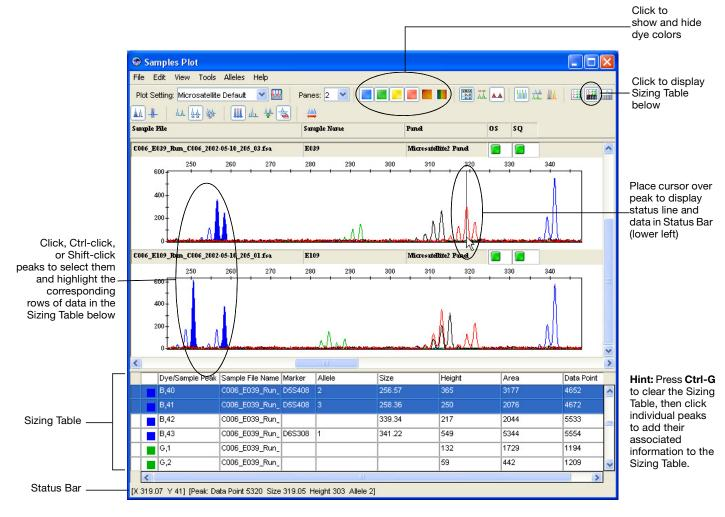


Figure 2-2 Examining and comparing data from different sample files in the Samples Plot



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 **Microsatellite Kit** you created on page 11.
- 3. Click | (Bins ▶ New Bin Set).
- **4.** In the New Bin Set dialog box, type **Microsatellite Bin Set** for the Bin Set Name, then click **OK**.



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

Adding Reference Data to a Panel and Bin Set

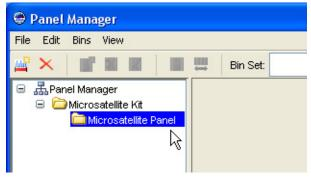
Note: You can add all or only a subset of the sample files in a project as reference data. Because there are a limited number of sample files, you will use all of them for reference data. If you have a large number of sample files, you could pick a subset of sample files that you believe to contain all of the alleles that exist in the sample files.

To add reference data to the Microsatellite Panel and Microsatellite Bin Set:

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



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

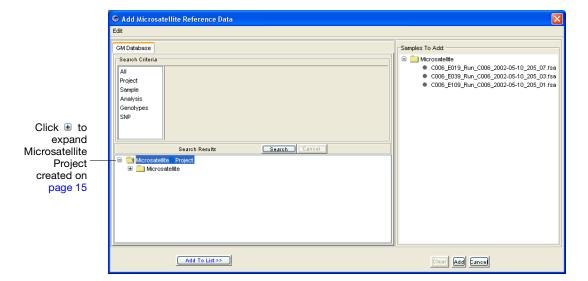


All the markers you created display on the right pane of the Panel Manager (Figure 2-3 on page 39).

3. Click **№** (Bins > Add Reference Data).

The Add Microsatellite Reference Data dialog box opens displaying the Microsatellite 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. Expand the **Microsatellite Project**, select the **Microsatellite** folder, click **Add to List**, then click **Add**.

All the sample files in the Microsatellite Project are added as reference samples to the Microsatellite Panel and appear in the lower half of the Navigation Pane in the Panel Manager (Figure 2-3 on page 39).

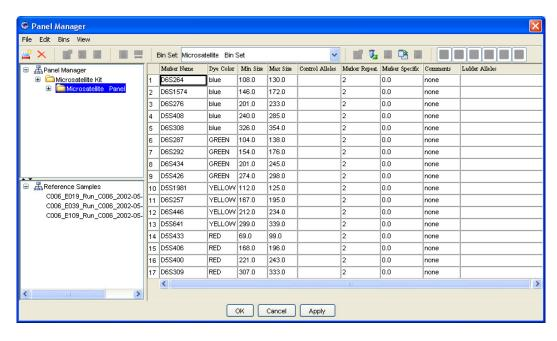
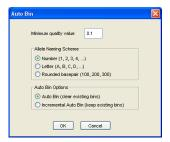


Figure 2-3 Panel Manager displaying markers and reference sample files for selected panel

Generating Bins using Auto Bin

To generate bins using the Auto Bin feature:

- 1. Make sure the Microsatellite Panel is selected in the Navigation Pane and Microsatellite Bin Set is selected in the Bin Set drop-down list, then click (Bins > Auto Bin).
- **2.** In the Auto Bin dialog box:
 - Leave Minimum quality value set to **0.1**
 - Select Number for Allele Naming Scheme
 - Select **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 Microsatellite 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*.

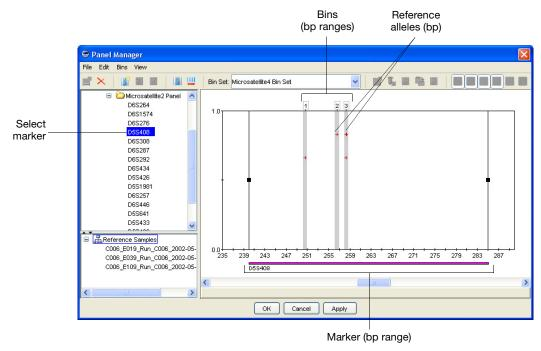


Figure 2-4 Selecting a marker

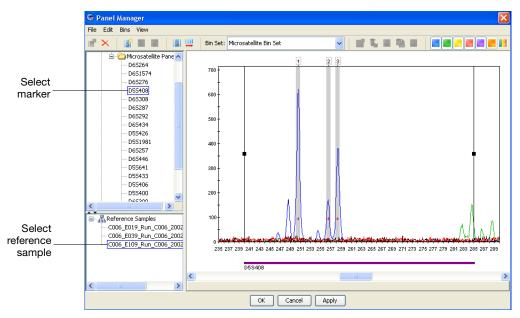


Figure 2-5 Selecting a marker and reference sample

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 Microsatellite Kit and Microsatellite 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

- 1. Select the marker in the upper Navigation Pane, then click (Bins > Add Bin).
- 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.

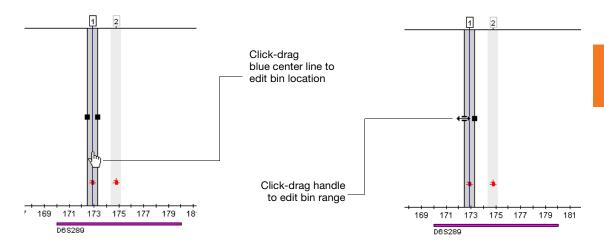


Figure 2-6 Editing a bin graphically

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 44).
- **2.** Click-drag the left or right handles that define the marker range.

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

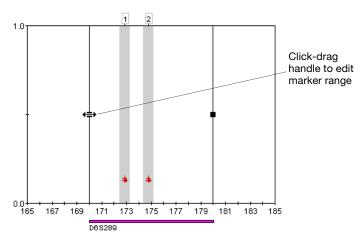


Figure 2-7 Editing a marker

Deleting a Marker From a Panel

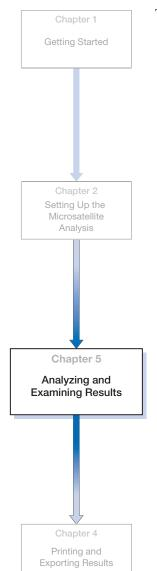
- 1. Select the marker in the upper Navigation Pane.
- 2. Click × (Edit → Clear Marker).

Next Steps

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

3

Analyzing and Examining Results



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

Editing the Analysis Method

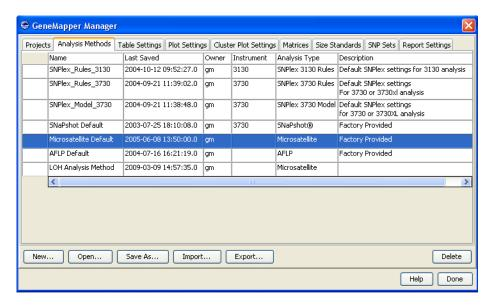
Overview

When you created the Microsatellite Analysis Method on page 18, 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® Software to make allele calls when you analyze the sample files in the project.

Editing the Analysis Method

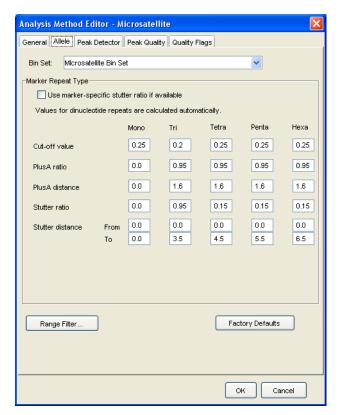
To edit the Microsatellite Analysis Method:

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



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

4. In the Analysis Method Editor, select the **Allele** tab, select **Microsatellite 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 18).

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[®] 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 Microsatellite 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

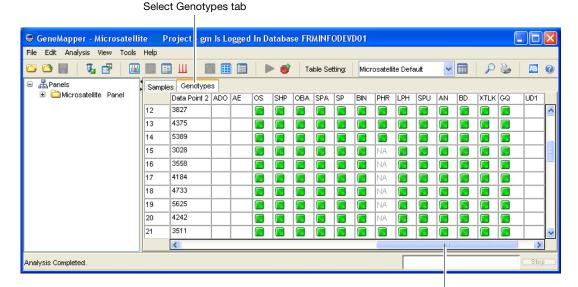
Overview

To examine the sizing and genotyping results:

- Review the SQ, associated PQVs, size standard, sample information, and samples plots (described on pages 28 through 34)
- 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 the Genotypes Plot window (page 52)
- View project alleles (page 52)

Reviewing the GO and POVs

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 .

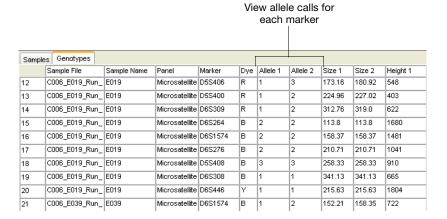
Investigating Yellow A and Red GQs

IMPORTANT! When analyzing your own data, you may find the GQ to be △ (Check) or ♠ (Low Quality) and the contributing PQVs (AN, BD, BIN, CC, 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.

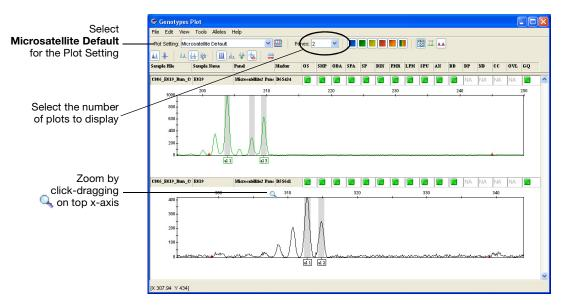


Viewing Genotype Plots

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 Zoom To , type range, then click OK .
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 Zoom To , type maximum, optionally, select Apply to all electropherograms , then click OK .
Unzoom	Double-click the x-axis or y-axis.
	or
	Right-click the x-axis or y-axis, then select Full View .

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*[®] *Software Online Help*.

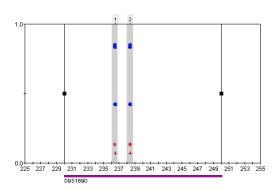
When done viewing the Genotypes Plot, click to close the window.

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 **Microsatellite Kit**, expand the **Microsatellite 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.





Printing and Exporting Results



This	chanter	includes:
1 1113	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 Analysis > Display 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*[®] *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 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	В
adding bins to markers 42 sample files to project 15 size standard labels 31 Advanced peak detection algorithm 21 allele calls, reviewing 50 Allele tab of Analysis Method Editor 20, 47 alleles, project 52 analysis method creating 18 editing 46 exporting 56 saving 24 selecting bin set 20, 47	bin adding to a marker 42 creating manually 36 definition 5, 36 deleting 43 editing 42, 43 generating using Auto Bin 40 reviewing 40 bin set accepting 42 creating 36 definition 5, 36 exporting 56 importing 36 selecting in analysis method 20, 47
Analysis Method Editor Allele tab 20, 47 General tab 19 Peak Detector tab 21 Peak Quality tab 22 Quality Flags tab 23 analysis parameters definition 5, 18 setting 18 analysis. See microsatellite analysis analyzing a project 27, 48 Applied Biosystems contacting viii customer feedback on documentation viii Information Development department viii Technical Support viii assumptions for using this guide v Auto Bin 40	C conventions bold text v for describing menu commands v IMPORTANTS! vi in this guide v italic text v Notes vi user attention words vi creating analysis method 18 bin set 36 bins 40 kit 11 markers 11 panel 11 project 15 reports 56 customer feedback, on Applied Biosystems

documents viii	G
D	General tab of Analysis Method Editor 19 genotype quality. See GQ
data See example data See reference data See sample files	Genotypes Plot window displaying 50 examining data in 52 exporting 55
deleting bins 43 markers 44 size standard labels 31	printing 54 zooming 51 Genotypes tab 49 Genotypes Table 34
documentation, related vii	GQ GQ
E	investigating 49 reviewing 49
editing analysis method 46 bins 42, 43 markers 44 size standard labels 31 example data	H help, online, accessing vii, viii heterozygous min peak height 22 hiding dye color peaks 34, 52
instrument used 2 marker information 3, 14 overview 2 size standard for 2	homozygous min peak height 22
exporting analysis methods 56 bin sets 56 Genotypes Plot window 55 kits 55 panels 55 plot settings 56	importing bin set 36 panels 11 Information Development department, contacting viii instruments
projects 56 report settings 56 reports 56 results 55	compatible with microsatellite analyses 2 for example data 2 italic text, when to use v
Samples Plot window 55 size standards 56 table settings 56	kits compatible with microsatellite analyses 2 creating 11
Files, See comple files	definition 5, 11 exporting 55
files. See sample files fill down 25	type 11

M	Editor 22
markers	plot setting 33, 51, 56
adding bins to 42	Plot window
creating 11	See Genotypes Plot window
definition 2, 5, 11	See Samples Plot window
deleting from a panel 44	plots. See sample plots
editing 44	primers, custom 2
for example data 3, 14	printing
plots, viewing 50 reviewing 40	Genotypes Plot window 54
viewing plots 50	reports 54
menu commands, conventions for	results 54
describing v	Samples Plot window 54
method. See analysis method	project
-	adding sample files to 15
microsatellite analysis	analyzing 27, 48
compatible instruments 2 compatible kits 2	creating 15 exporting 56
definition 2	saving 26
flowchart 4	setting analysis parameters 18
markers 2	setting table settings 18
setting up 9	project alleles
min peak height ratio 22	viewing 52
MSDSs, obtaining viii	
	Q
0	Quality Flags tab of Analysis Method
online help, accessing vii, viii	Editor 23
	Editor 25
overriding SQ 31	R
5	n
P	reference data
Panel Manager	adding to a panel 37
opening 11	icon 48
panels	report settings, exporting 56
adding reference data 37	reports
creating 11	creating 56
definition 5, 11	exporting 56
exporting 55 importing 11	printing 54
selecting for analysis 24	results
peak detection algorithm 21	exporting 55 printing 54
Peak Detector tab of Analysis Method	printing 34
Editor 21	S
Peak Quality tab of Analysis Method	sample files

adding to project 15	investigating 28
location 2	overriding 31
plots, viewing 32, 50	reviewing 28
sorting 28, 49	score 29
viewing information on 32	Status Bar 27
viewing plots 32, 50 zooming 33, 51	
	T
sample plots	table settings 18, 25, 56
viewing 32, 50	
zooming 33, 51	Technical Support, contacting viii
samples sorting 28, 49	text conventions v
<u> </u>	training, information on viii
zooming 33, 51	troubleshooting the size standard 31
Samples Plot window	S
displaying 32	U
examining data in 34 exporting 55	O .
printing 54	unzooming 33, 51
zooming 33	user attention words, described vi
Samples tab 18	
_	X
saving	
analysis method 24 project 26	x-axis
	scale 34, 52
setting up analysis 9	zooming 33, 51
showing and hiding dye color peaks 34, 52	
size calling curve 30	Y
Size Match Editor 29	y-axis
size quality. See SQ	scale 34, 52
size standard	zooming 33, 51
custom 31	,
examining 29	Z
exporting 56	-
for example data 2	zooming
labels, adding 31	Genotypes Plot window 51
labels, deleting 31	Samples Plot window 33
labels, editing 31	unzooming 33, 51
selecting for analysis 24	x-axis 33, 51
troubleshooting 31	y-axis 33, 51
Sizing Table 34	
software	
starting and logging in 5	
terms defined 5	
sorting samples 28, 49	
SQ	

