

GeneScan Analysis Software

Version 3.1

User's Manual

© Copyright 1998, The Perkin-Elmer Corporation

This product is for research purposes only.

GeneScan, Genotyper, and Perkin-Elmer are registered trademarks of The Perkin-Elmer Corporation.

BioLIMS is a trademark of Molecular Informatics, Inc.

ABI, ABI PRISM and the ABI PRISM design, Applied Biosystems, GenoPedigree, PE, and PE Applied Biosystems are trademarks of The Perkin-Elmer Corporation.

Sybase is a registered trademark and Sybase SQL Server are trademarks of Sybase, Inc.

Oracle is a trademark of Oracle Corporation.

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

Contents

1	<i>GeneScan Analysis Software Overview</i>	1-1
	Introduction	1-1
	In This Chapter	1-1
	What's New in the GeneScan Analysis Software	1-2
	BioLIMS Support	1-2
	Fifth Dye	1-2
	Works with 96-Lane Gels	1-2
	Displaying Off-Scale Data	1-3
	Manual Tracker Interface Changes	1-3
	Printing from Sample File Window	1-3
	New Neural Net Tracker	1-3
	Lane Extraction	1-4
	About This Manual and Other Instruments	1-5
	ABI 373	1-5
	ABI 373 XL Upgrade Users	1-5
	96-lane Upgrade Kit	1-5
	Manual Text	1-5
	User Attention Words	1-6
	Using the Macintosh Computer	1-7
	Macintosh Computer Vocabulary and Operations	1-7
	Guidelines for Optimal Performance	1-7
	Macintosh Computer Terms Used in This Manual	1-8
	Registering the Software	1-11
	License and Warranty	1-11
	Registering Your Software	1-11
	Installing the Software	1-12
	System Specifications	1-12

Turning On Virtual Memory	1-13
RAM Required to Run the Software	1-14
Installing the Software	1-14
Program Files Installed Diagram	1-15
Contents of Folders	1-15
GeneScan Analysis Software Files	1-16
Table of Files	1-16
Sample File Database Support	1-19
Technical Support	1-20
To Reach Us On the Web	1-20
Hours for Telephone Technical Support	1-20
To Reach Us by Telephone or Fax in North America	1-20
Fax-on-Demand	1-22
To Reach Us by E-Mail	1-23
Regional Offices Sales and Service	1-23

2 How to Process and Edit the Gel File 2-1

Introduction	2-1
In This Chapter	2-1
Topics in This Chapter	2-1
About Gel Files	2-2
Instruments Producing Gel File	2-2
96-Lane Gels Capability	2-2
Gel File Contents	2-2
How GeneScan Tracks a Gel File	2-2
How to Set Gel Processing Preferences	2-3
Introduction	2-3
Displaying the Gel Preferences Dialog Box	2-3
Auto-Launch Processing	2-4
Image Generation Defaults	2-4
Lane Extraction	2-7
Comb Type	2-10
How to Display the Gel File	2-11
Introduction	2-11

Displaying the Gel File	2-11
About the Gel File Window	2-12
Differences Between Gel File Windows	2-12
Gel File Window Diagram	2-13
Gel File Window Buttons	2-14
Gel File Window Described	2-15
How to Adjust the Gel Image	2-18
Introduction	2-18
Displaying/Hiding a Dye Color	2-18
Changing Dye Indicators	2-19
Adjusting the Contrast	2-20
Removing Gel Contrast Changes	2-22
Regenerating the Gel Image	2-23
Installing New Matrix Information	2-24
Verifying Gel Information	2-26
Where the Information is Archived	2-26
Data Collection Run Error Log	2-26
Gel Info Window	2-26
Gel Image	2-27
Sample Sheet	2-28
Displaying the Sample Sheet	2-29
What is the Sample Sheet	2-29
Using the Copy of the Sample Sheet	2-29
Displaying the Copied Sample Sheet	2-29
Displaying a New Sample Sheet	2-29
About the Sample Sheet	2-31
Sample Sheet Example	2-31
Sample Sheet Described	2-31
Working with the Sample Sheet	2-34
Editing Collection Information	2-34
Changing Column Widths	2-35
Editing the Sample Sheet	2-35
Printing the Sample Sheet	2-36
Installing a New Sample Sheet	2-36
Importing Data to Sample Sheets	2-37

Exporting Data From Sample Sheets	2-37
About Displaying Regions of Off-Scale Data.	2-38
Why Displaying Off-Scale Data is Important	2-38
Off-Scale Detection and GeneScan	2-38
Displaying Off-Scale Data	2-38
Gel File Window Example	2-40
Electropherogram Examples	2-41
Working with Lane Markers	2-43
Introduction	2-43
Topics in This Section	2-43
Moving Between Lanes and Channels	2-44
Using Control Points	2-45
Adding and Deleting Control Rows	2-46
What Marking Lanes as Used/Unused Means	2-47
Marking Lanes as Used or Unused	2-47
Moving Misplaced Lane Markers	2-48
Marking and Unmarking Lanes for Extraction.	2-50
Lane Marker Rules	2-50
About Marking Lanes for Extraction	2-50
Marking All Lanes for Extraction	2-51
Unmarking All Lanes for Extraction	2-51
Marking a Single Lane for Extraction	2-52
Unmarking a Single Lane for Extraction	2-52
Working with Tracker Lines	2-53
Topics in This Section	2-53
Hiding/Showing Tracker Lines	2-53
Why Positioning Each Lane is Important	2-54
Causes to Misinterpreting Lane Positions	2-54
Verifying Optimal Channel Tracking	2-55
Reviewing Tracker Lines	2-55
Moving Tracker Lines	2-56
Deselecting a Lane	2-56
Interpolating Tracker Lines.	2-57
Why Use Interpolation Mode.	2-57
Procedure	2-57

Tracking Lanes and Extracting Data	2-59
Ways to Track and Extract Data	2-59
If the Run Folder is Not Specified	2-59
Tracking the Gel Without Extracting Data	2-60
Tracking and Extracting Data	2-61
Extracting Data Without Changing the Current Tracker.	2-64
How the GeneScan Analysis Software Names Sample Files	2-67
Naming Process	2-67
Saving Gel Files After Editing Tracking	2-68
Introduction.	2-68
Saving Files After Editing Tracking	2-68

3 Analyzing Sample Files 3-1

Introduction	3-1
In This Chapter	3-1
About Sample Files	3-2
What Sample Files Contain	3-2
How GeneScan Generates Sample Files.	3-2
Ways to Generate Sample Files	3-3
How GeneScan Analyzes Sample Files	3-3
Opening Sample Files	3-4
Introduction.	3-4
Procedure	3-4
About The Sample File Window	3-5
What it Displays	3-5
Five Views.	3-5
Sample Results View	3-6
What it Displays	3-6
Displaying the View	3-6
Sample Results View Example	3-6
Description of Columns	3-7
Differences From the Results Display	3-7
Sample Info View.	3-8
What it Displays	3-8

Displaying the View	3-8
Sample Info View Example	3-9
Description of Information.	3-9
Size Curve View	3-13
What it Displays.	3-13
Displaying the View	3-13
What the Size Curve Displays	3-14
Curves Described.	3-14
Raw Data View	3-15
What it Displays.	3-15
Displaying the View	3-15
Raw Data Example.	3-16
What to Evaluate	3-16
EPT Data View	3-17
What it Displays.	3-17
Displaying the View	3-17
EPT Data Example.	3-18
Colored Lines Described	3-18
Analyzing a Sample File.	3-19
Introduction	3-19
Procedure	3-19
Installing a New Matrix File	3-20

4 *Creating a Project*4-1

Introduction.	4-1
In This Chapter.	4-1
Automatic Analysis and Project Creation Process	4-2
Introduction	4-2
Process Using the ABI PRISM 310	4-2
Process Using the ABI 373 and the ABI PRISM 377	4-3
Setting Up for Automatic Analysis	4-4
Introduction	4-4
Procedure	4-4
Using a Project to Manage Sample Files	4-8

What is a Project	4-8
Why Create a Project	4-8
When GeneScan Creates Projects	4-8
Where to Store Projects	4-9
Working with Project Files	4-10
Opening an Existing Project	4-10
Creating a New Project	4-10
Unlocking Sample Files	4-12
Removing Samples from a Project	4-13
Finding Missing Sample Files	4-14
When are Files Considered Lost	4-14
When an Alert Appears	4-14
Searching for Missing Sample Files	4-14
Re-Establishing Links	4-15

5 Analyzing Project Files 5-1

Introduction	5-1
In This Chapter	5-1
Analyzing Project Files: About the Analysis Control Window	5-2
Introduction	5-2
Analysis Control Window Example	5-2
Analysis Control Window Description	5-3
Customizing the Display	5-4
Using the Analysis Control Window	5-4
Analyzing Sample Files: Using the Analysis Control Window	5-6
Introduction	5-6
Accessing Sample Files	5-6
Analyzing Sample Files	5-7
Specifying the Format for Printed Results	5-9
Displaying Size Standards and Analysis Parameters	5-11
Displaying Sample and Dye Information	5-12
Setting Sample File Sort Order	5-14
Setting Dye Indicator Preferences	5-15
About the Analysis Parameters	5-17

Introduction	5-17
Parameters Used During Automatic Analysis	5-17
Specifying Where Files are Stored.	5-17
Defining Analysis Parameters	5-18
Six Analysis Parameters.	5-18
Analysis Range Parameter Options	5-19
Data Processing Parameter Options	5-19
Peak Detection Parameter Options.	5-20
Size Call Range Parameter Options	5-21
Size Calling Method Parameter Options	5-21
Split Peak Correction Parameter Options	5-22
Using Analysis Parameter Files	5-24
In This Section	5-24
Analyzing Samples Using the Same Analysis Parameters.	5-24
Selecting Different Parameters for Analysis Samples	5-25
Displaying Default Parameters	5-26
Creating Custom Analysis Parameter Files	5-26
Changing an Existing Analysis Parameters File	5-27
Deleting Custom Analysis Parameters.	5-28
About Size Standards	5-29
What are Size Standards	5-29
Advantages of Using Size Standard.	5-29
Size Standards Provided.	5-29
When to Define Size Standards	5-29
If Split Peaks Appear	5-29
Specifying Where Files are Stored.	5-30
Defining the Size Standard	5-31
Two Ways to Define the Size Standard	5-31
Using the New Command	5-31
Using the Analysis Control Window	5-34
Using Size Standards	5-36
In This Section	5-36
Changing the Number of Peaks Detected	5-36
Editing the Size Standard Definition	5-36
Editing an Existing Standard	5-37

Deleting an Existing Standard	5-39
Analyzing Samples Using the Same Standard	5-39
Selecting Separate Standards for Samples	5-40
Defining Folder Locations	5-41
Introduction	5-41
Storing Matrix Files	5-41
Specifying File Locations	5-42

6 *Making a Matrix File* 6-1

Introduction	6-1
In This Chapter	6-1
About Matrix Files	6-2
Introduction	6-2
Matrix File Definition	6-2
Multicomponent Definition	6-2
Why is a Matrix File Necessary	6-3
When to Create a Matrix File	6-3
Sample Files Using Matrix File	6-4
Applying a Matrix File	6-4
When to Assign a Matrix File	6-4
Limitations to Matrix Files	6-5
When to Create New a Matrix File	6-5
Considerations Before Making a Matrix File	6-6
Where to Store Matrix Files	6-6
Process of Creating a New Matrix File	6-7
Process Diagram	6-7
Steps to Creating a New Matrix	6-8
The Dye Matrix Standard Kits	6-8
Table of Kits	6-8
Preparing Matrix Standards for the ABI PRISM 310	6-9
About Formamide and Samples in Formamide	6-9
Deionizing Formamide	6-9
Preparing the Formamide-Size Standard Mix	6-10
Preparing Matrix Standard Samples	6-10

.....	6-11
Preparing Matrix Standards for the ABI 373 and ABI PRISM 377	6-12
Introduction	6-12
Procedure	6-12
Table of Matrix Standards	6-12
Loading and Running Dye Standards for the ABI PRISM 310	6-13
Introduction	6-13
Creating a GeneScan Sample Sheet	6-13
Creating a GeneScan Injection List	6-14
Starting the Data Collection Software	6-15
If the Run was Cancelled	6-15
Run Time	6-16
Assigning Matrix File	6-16
Loading and Running Dye Standards for the ABI PRISM 377	6-17
Introduction	6-17
Creating a GeneScan Sample Sheet	6-17
Loading Matrix Standards	6-18
Running the Matrix Standards	6-19
Loading and Running Matrix Samples on the ABI 373 (non XL)	6-20
Introduction	6-20
Loading the Matrix Standards	6-20
Completing the Sample Sheet in the 672 Software	6-21
Running the Gel	6-22
Completing the GeneScan Sample Sheet	6-23
Loading and Running Matrix Standards on the ABI 373 with XL Upgrade	6-24
Introduction	6-24
Completing the GeneScan Sample Sheet	6-24
Loading the Matrix Standards	6-25
Running the Matrix Standard	6-26
Generating Matrix Sample Files for the ABI 373 and ABI PRISM 377 ...	6-27
Who Should Use This Step	6-27
Verify Tracking	6-27
Generating Sample Files	6-27
Choosing a Scan Range for the Matrix Calculation	6-29
Introduction	6-29

Viewing the Raw Data	6-29
What to Look For in the Raw Data Display	6-30
Choosing a Scan Range	6-30
Eliminating Primer Peaks	6-31
Generating a New Matrix File	6-32
Procedure	6-32
Matrix File Example	6-33
Saving and Naming the Matrix File.	6-34
Introduction.	6-34
Naming Considerations.	6-34
Saving the Matrix File	6-34
Where to Store the Matrix File	6-34
When to Make Copies of a Matrix	6-34
Assigning the Matrix File to Sample Files	6-35
Introduction.	6-35
Procedure	6-35
Evaluating the Matrix File.	6-37
Introduction.	6-37
Procedure	6-37
Causes for Bad Matrix Files	6-38
If an Error Message Appears.	6-38
Two Causes of Bad Matrix Files.	6-38

7 Evaluating Analysis Results. 7-1

Introduction	7-1
In This Chapter	7-1
Process of Evaluating Analysis Results	7-2
Steps to Evaluating Analysis Results	7-2
Ways to Display Analysis Results	7-3
Two Ways to Display Analysis Results	7-3
The Results Display Window	7-4
Introduction.	7-4
Displaying the Window.	7-4
Results Control Window Description	7-5

Using the Results Control Window	7-7
Selecting Display Format	7-7
Selecting Electropherogram Panels	7-7
Selecting Samples to Display	7-8
Creating Tiled Electropherogram Displays	7-10
Deselecting Samples	7-11
Removing Samples	7-12
Displaying the Results	7-12
Printing the Results	7-12
Changing How the Results are Displayed and Printed	7-13
Procedure	7-13
The Sample Results View	7-16
About the View	7-16
For More Information	7-16
Updating the Results	7-16
Re-Analyzing the Data	7-16
Saving and Renaming the Results Control Format	7-17
Introduction	7-17
Important Considerations	7-17
Saving the Display Format	7-17
Working with a Previously Saved Displays	7-18
Renaming the Current Results Display	7-19
About Electropherogram and Tabular Data Displays	7-20
Introduction	7-20
How the Window is Divided	7-20
What Tabular Data Contains	7-20
How Electropherogram Panels are Sized	7-20
Peaks Visible in Electropherogram Not Tabular Data	7-21
For More Information	7-21
Displaying Electropherogram and Tabular Data	7-22
Displaying Electropherogram and Tabular Data	7-22
Tabular Data and Electropherogram Example	7-22
Table Describing Columns	7-23
Highlighting Information	7-23
Adjusting Window Size	7-24

Hiding Selected Rows of Data	7-24
Limited the Rows Displayed.	7-24
Displaying Electropherogram Data	7-25
Definition	7-25
Base Pairs Versus Data Points.	7-25
Procedure	7-25
Electropherogram Example	7-26
Descriptions of the Electropherogram	7-26
Working with Electropherogram Data.	7-27
In This Section	7-27
Displaying X-and Y-Axis Positions	7-27
Moving the Electropherogram	7-28
Changing the Dye Color	7-28
Displaying Peak Positions.	7-29
Highlighting Peaks	7-29
Using Legends to Change the Display	7-30
Scrolling the Display.	7-30
Zooming In and Out	7-31
Showing Off-Scale Data	7-32
Electropherogram Displaying Off-Scale Data	7-34
Electropherogram Displaying the Flat Topped Effect	7-35
Showing Data by Fragment Size.	7-35
Changing the Horizontal Scale	7-37
Changing the Vertical Scale	7-38
Assigning Standard or Custom Colors	7-39
How to Define Custom Colors.	7-40
Introduction.	7-40
Why Change Colors in the Electropherogram	7-40
Saving the Display Format	7-40
Defining Custom Color for All Electropherogram	7-41
Defining Individual Plot Colors	7-42
How to Change the Dye Scale	7-44
What the Dye Scale Defines	7-44
Increasing the Dye Scale Example	7-44
Changing the Dye Scale of an Electropherogram.	7-44

Changing the Dye Scale Preferences	7-45
Process of Verifying Results	7-46
Introduction	7-46
Steps to Verifying Size Calculation	7-46
How to Verify Size Calculations	7-48
Introduction	7-48
Verifying for the GeneScan-350 Standard	7-48
Evaluating for Multiple Internal Size Standards	7-50
Using the Analysis Log	7-51
What is the Analysis Log	7-51
Displaying the Analysis Log	7-51
What to Evaluate	7-52
Removing Information from the Analysis Log	7-52
Closing the Analysis Log	7-52
How to Verify Peak Detection	7-53
Introduction	7-53
Verifying Peak Detection	7-53

8 *Saving, Archiving and Copying Files8-1*

Introduction	8-1
In This Section	8-1
Saving GeneScan File	8-2
Why Save	8-2
Saving Projects	8-3
Saving Sample Files	8-3
Saving Gel Files	8-4
Saving Results Displays	8-5
How to Archive Sample and Gel Files	8-6
Introduction	8-6
Archiving Sample and Gel Files	8-6
Transferring Data to Other Applications	8-7
Other Applications	8-7
Cutting and Pasting Tabular Data	8-7
Creating a Text File	8-8

9	<i>Printing Results</i>	9-1
	Introduction	9-1
	In this Chapter	9-1
	About Printing	9-2
	Ways You Can Print	9-2
	If You Get Unexpected Results	9-2
	Automatically Printing Run Results	9-3
	Introduction	9-3
	From the Data Collection Software	9-3
	From the GeneScan Analysis Software	9-3
	Printing the Gel Image	9-4
	About Printing the Gel Image	9-4
	Procedure	9-4
	Printing Selected Sample Files	9-5
	Introduction	9-5
	Setting Printing Options	9-5
	From the Results Control Window	9-5
	From the File Menu	9-6
	Printing Supporting Information	9-6
	Introduction	9-6
	Procedure	9-6
A	<i>Using GeneScan with the ABI 373</i>	A-1
	Introduction	A-1
	In This Appendix	A-1
	Processing Using the ABI 373	A-2
	Introduction	A-2
	ABI 373 Versus ABI PRISM 377 Data	A-2
	Using the GeneScan Analysis Software Before Version 2.0	A-2
	Using the GeneScan Analysis Software Version 2.0 or Later	A-2
	Processing Steps	A-3
	Comparison of GeneScan Analysis	A-3
	Setting Gel Processing Parameters	A-5

Introduction	A-5
Procedure	A-5
Opening the Collection File	A-7
Introduction	A-7
Procedure	A-7
Completing the Sample Sheet.	A-9
Introduction	A-9
Procedure	A-9
Tracking Lanes	A-11
Introduction	A-11
Process	A-11
Procedure	A-11
Generating Sample Files.	A-12
When to Generate Sample Files.	A-12

B GeneScan Size Standards B-1

Introduction.	B-1
In This Appendix	B-1
GeneScan 350 Size Standard	B-2
What To Use It For.	B-2
How It Is Prepared	B-2
GeneScan 350 Molecular Lengths	B-2
Running Under Denaturing Conditions	B-2
Double-Stranded GeneScan 500 Fragments	B-3
GeneScan-400HD Size Standard	B-4
What To Use It For.	B-4
Special Uses	B-4
How It Is Prepared	B-4
Fragment Lengths	B-4
Denaturing Electropherogram	B-5
Non-denaturing Electropherogram.	B-5
GeneScan 500 Size Standard	B-6
What To It Use For.	B-6
How It Is Prepared	B-6

GeneScan 500 Molecular Lengths	B-6
Running Under Denaturing Conditions	B-6
Double-Stranded GeneScan 500 Fragments	B-7
GeneScan 1000 Standard	B-8
How It Is Prepared	B-8
GeneScan 1000 Molecular Lengths	B-8
Running Under Native Conditions	B-9
GeneScan 2500 Standard	B-10
How It Is Prepared	B-10
What To Use It For	B-10
Running Under Native Conditions	B-10
GeneScan 2500 Molecular Lengths	B-11

C Size Calling Methods C-1

Introduction	C-1
In This Appendix	C-1
Least Square Method	C-2
About This Method	C-2
Advantages	C-2
Cubic Spline Interpolation Method	C-4
About This Method	C-4
Possible Local Sizing Inaccuracy	C-4
Local Southern Method	C-5
About This Method	C-5
The Equation	C-5
How This Method Works	C-6
Global Southern Method	C-7
About This Method	C-7
The Equations	C-7
How This Method Works	C-8

D Troubleshooting the GeneScan Software D-1

Introduction	D-1
------------------------	-----

In This Appendix	D-1
Troubleshooting Projects and Results	D-2
Table Description	D-2
Troubleshooting Gel Data	D-5
Table Description	D-5
Troubleshooting Genotyping Results	D-8
Table Description	D-8
GeneScan Error Messages	D-9
Introduction	D-9
GeneScan Analysis Software Crashes with BioLIMS	D-9
Analysis Log Error Messages	D-10
Error Messages When Defining Size Standards	D-11

E Using the BioLIMS Database E-1

Introduction	E-1
In This Appendix	E-1
About the GeneScan Analysis Software and the BioLIMS Database	E-2
What is the BioLIMS Database?	E-2
Accessing the BioLIMS Database	E-2
Before Using the BioLIMS Database	E-2
Modes in GeneScan Analysis Software	E-2
Comparing Modes	E-3
Configuring the BioLIMS Database Server	E-4
Sybase or Oracle?	E-4
Configuring for Sybase SQL Server Connection	E-4
Configuring for Oracle Server Connection	E-7
Switching Between Sample File and BioLIMS Mode	E-10
Introduction	E-10
Switching Modes	E-10
How to Access the BioLIMS Database	E-12
Introduction	E-12
Before Accessing a Sybase Database	E-12
Accessing the BioLIMS Database	E-13
About Server Names	E-17

Sybase or Oracle?	E-17
How Names are Recognized.	E-17
Sybase SQL Server Examples	E-17
Oracle Server Examples	E-18
Using the Collection Browser Window	E-20
Applications that Use the Collection Browser Window	E-20
Ways to Search the Database	E-20
In This Section	E-20
Displaying the Collection Browser Window.	E-21
Collection Browser Window Example	E-22
Parts of the Collection Browser Window	E-22
Collection Search Criteria.	E-24
Fragment Search Criteria	E-25
Searching the BioLIMS Database.	E-28

F Troubleshooting the BioLIMS Database. F-1

Introduction	F-1
In This Appendix.	F-1
If the BioLIMS™ Preference Page Does Not Appear	F-2
Problem.	F-2
Solution.	F-2
About Troubleshooting the Client to Sybase Connection	F-3
Introduction.	F-3
SybPing and Telnet	F-3
Troubleshooting Process	F-4
Troubleshooting Flow Chart.	F-4
Procedures for Troubleshooting the Client to Sybase Connection	F-6
Introduction.	F-6
Troubleshooting from the Macintosh Client.	F-6
Required Sybase Extension Files	F-11
Required Oracle Extension Files	F-11
Troubleshooting from the Unix (Sybase) Server	F-12
About Troubleshooting the Client to Oracle Connection	F-14
Introduction.	F-14

Telnet	F-14
Troubleshooting Process	F-15
Troubleshooting Flow Chart	F-15
Procedures for Troubleshooting the Client to Oracle Connection	F-17
Introduction	F-17
Troubleshooting from the Macintosh Client	F-17
Troubleshooting from the Unix (Oracle) Server	F-22

G License and Warranty G-1

Software License and Warranty	G-1
Perkin-Elmer Software License and Limited Product Warranty	G-1
Copyright	G-1
License	G-1
Restrictions.	G-2
Limited Warranty	G-2
Limitation Of Liability.	G-2
Term	G-3
Miscellaneous.	G-3

Glossary

Index

GeneScan Analysis Software Overview

1

Introduction

In This Chapter This chapter provides a general introduction to the ABI PRISM® GeneScan® Analysis Software. It provides information about the organization of this manual and instructions on how to get help from PE Applied Biosystems.

Topics in this chapter include the following:

Topic	See page
What's New in the GeneScan Analysis Software	1-2
About This Manual and Other Instruments	1-5
Using the Macintosh Computer	1-7
Registering the Software	1-11
Installing the Software	1-12
GeneScan Analysis Software Files	1-16
Technical Support	1-20

Note Already familiar with previous versions of the GeneScan Analysis Software and want to know what is new and different in this version? Turn to "What's New in the GeneScan Analysis Software" on page 1-2.

What's New in the GeneScan Analysis Software

BioLIMS Support GeneScan Analysis Software version 3.1 supports the BioLIMS™ Genetic Information Management System v. 2.0. In the BioLIMS mode, the GeneScan Analysis Software extracts sample data from gel files and writes it to the database. The contents of the gel file are not saved to the database. The GeneScan Analysis Software can read and write sequence-fragment to the database just as it writes sample files to the Macintosh® hard disks.

Both Oracle® and Sybase® databases are supported.

For information about using the GeneScan Analysis Software in BioLIMS mode, see “Using the BioLIMS Database” on page E-1.

Fifth Dye The GeneScan Analysis Software version 3.1 can display and analyze raw data from up to five dye colors. This feature has been implemented in the software to support future developments in the dye chemistries.

The current maximum number of dyes available is four. The availability of a fifth dye for the GeneScan Analysis Software is dependent on the progress made in the dye development chemistry and can not be accurately projected at this point. You can receive more information on the availability of additional dyes for the GeneScan Analysis Software when they become available through our web site (see “To Reach Us On the Web” on page 1-20).

Works with 96-Lane Gels The GeneScan Analysis Software version 3.1 can open and analyze 96-lane gels. For more information, refer to the *ABI PRISM 377 DNA Sequencer 96-Lane Upgrade User's Manual* (P/N 4305423).

Displaying Off-Scale Data

A new feature has been added in version 3.1 to display offscale regions. By clicking the OS button on the tool bar when the gel image is displayed, if the gel contains offscale data, then this data appears as white bands on the gel image. Off-scale data can also be viewed while in Results Display by selecting Show Offscale Regions from the View menu.

Note To see off-scale portions in the gel image, click the Vertical Expand button. The resolution of the full display is low and does not show the off-scale data.

Manual Tracker Interface Changes

In version 3.1, the new tracker lines consist of control points. The new interface allows the selection and movement of multiple control points simultaneously, interpolation of tracker line shapes between two lanes, multiple lane selection, and movement.

Printing from Sample File Window

In version 2.1, only the electropherogram display is printed when printing from the Sample File window.

The following displays can now be printed also:

- ◆ Sample Info View
 - ◆ Size Curve View
 - ◆ Raw Data View
 - ◆ EPT Data View
-
-

New Neural Net Tracker

About the Neural Net Tracker

This was the most significant improvement in the version 3.1 release of GeneScan Analysis Software.

- ◆ The Neural Net Tracker program uses a neural net-based algorithm to automatically track gel lanes. The Neural Net Tracker has been taught how to recognize bands and how to track curved lanes.

IMPORTANT Before auto-tracking the first time, you need to verify that the comb type is set correctly. The default comb type is Square.

- ◆ The Neural Net Tracker program exists as a separate program within the GeneScan Analysis Software folder. Also associated with the Neural Net Tracker program are a set of Tracker settings files that have been optimized for number of lanes and comb-types.

- ◆ The Neural Net Tracker program is “headless.” This means that although it stands as a separate program file, it does not have a user interface. The Tracker program is opened automatically from within the GeneScan Analysis Software.

IMPORTANT The gel file must be multicomponented using the correct matrix in order to be auto-tracked.

Table of Tracking Times

Tracking times depend upon the number of lanes, channels, and scans in the gel file. Consult the table below to estimate gel tracking times for your GeneScan Analysis Software.

Number of Lanes	Number of Channels	Number of Scans	Size of Gel (MB)	Time (min) for CPU/Speed			
				7200/90 ^a	4400/200 ^a	9500/200 ^b	G3/266 ^a
36	194	9152	21.6	15	9	c	5
36	194	11424	26.8	20	11	c	6
48	388	4516	21.5	26	10	c	5
48	388	7768	36.1	38	17	c	8
64	388	4565	21.7	26	10	c	6
64	388	5708	26.8	30	13	c	6
96	480	6840	39.5	46	24	c	13

a. 7200, 4400, G3:32MB + 10MB VM

b. 9500: 64MB + 10MB VM

c. Could not test this system

Lane Extraction In version 3.1, you can extract lanes using weighted averaging or pre-averaging offscale detection. These options are available in the Lane Extraction section of the Gel Preferences dialog box.

For more information, see “Lane Extraction” on page 2-7.

About This Manual and Other Instruments

ABI 373 GeneScan Analysis Software works version 3.1 with ABI™ 373 data in the same way as the it works with ABI PRISM® 377 data.

If you use the GeneScan® 672 Software to collect data on the ABI 373, you must transfer your data to a Power Macintosh® computer, and perform certain extra steps. These extra steps are described in Appendix A, "Using GeneScan with the ABI 373."

ABI 373 XL Upgrade Users The ABI™ 373 DNA Sequencer with XL upgrade updates the data collection interface to match the ABI PRISM 377, increases the maximum image resolution to 388 channels, and increases the maximum number of lanes to 66.

Note ABI 373 with XL upgrade users do not need to follow the procedures described in this Appendix A, "Using GeneScan with the ABI 373."

96-lane Upgrade Kit The ABI PRISM® 377 DNA Sequencer 96-lane upgrade kit enhances both the capabilities of the 377 DNA sequencer to support up to 96 lanes for using the GeneScan Analysis Software.

Manual Text **Manual Text Applies to other ABI PRISM Genetic Analysis Instruments**

The GeneScan Analysis Software and this manual apply to three different genetic analysis instruments supported by PE Applied Biosystems.

Instrument	Description
ABI PRISM® 310 Genetic Analyzer	Analyzes one sample at a time using capillary electrophoresis technology. This instrument provides automatic sample loading while using a minimal amount of sample.
ABI 373 DNA Sequencer	Performs slab gel electrophoresis, allowing the user to analyze multiple samples on a gel.
ABI PRISM 377 DNA Sequencer	A high-throughput slab gel electrophoresis instrument, created to meet the needs of high-volume DNA sequencing or genetic analysis laboratories. Throughput is up to than four times that of the ABI 373.

**User Attention
Words**

The text of this manual includes the following user attention words to draw your attention to specific details of the information presented in this manual.

User Attention word	Description
Note	Used to call attention to information.
IMPORTANT	Indicates information that is necessary for proper operation of the software.

Using the Macintosh Computer

Macintosh Computer Vocabulary and Operations

To use the GeneScan Analysis Software, you should be familiar with the following basic Macintosh® computer vocabulary and operations:

Vocabulary and operations	Description
Using the mouse	Clicking and double-clicking, selecting and dragging.
Choosing commands	Using pull-down and pop-up menus, dialog boxes, radio buttons and checkboxes.
Working with windows	Opening and closing, resizing and repositioning, scrolling, understanding the active window.
Using the Macintosh computer hierarchical file system	Finding files and creating folders.

If you do not understand some or all of these, refer to the *Macintosh® System Software User Guide* or the *Owner's Guide* for more information.

Guidelines for Optimal Performance

Computers require regular attention and maintenance to operate efficiently and consistently. The ABI PRISM® software works with large files and accesses the hard disk often, so it is especially important that you follow maintenance procedures to minimize the occurrence of errors during operation.

Follow these general guidelines for optimal performance:

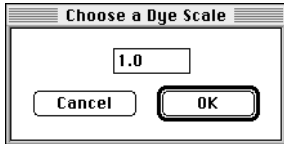
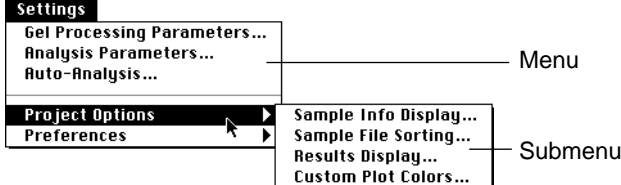
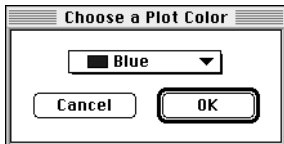
- ◆ Install only one Macintosh® computer system per hard disk.
- ◆ Back up all programs and files regularly.
- ◆ Remove nonessential files from the hard disk.
- ◆ Use a hard disk maintenance program regularly.
- ◆ Use discretion when adding software programs, especially Control Panel (CDEV) and Extension (INIT) files.

Refer to the *ABI PRISM Instrument User Manual* for more detail.

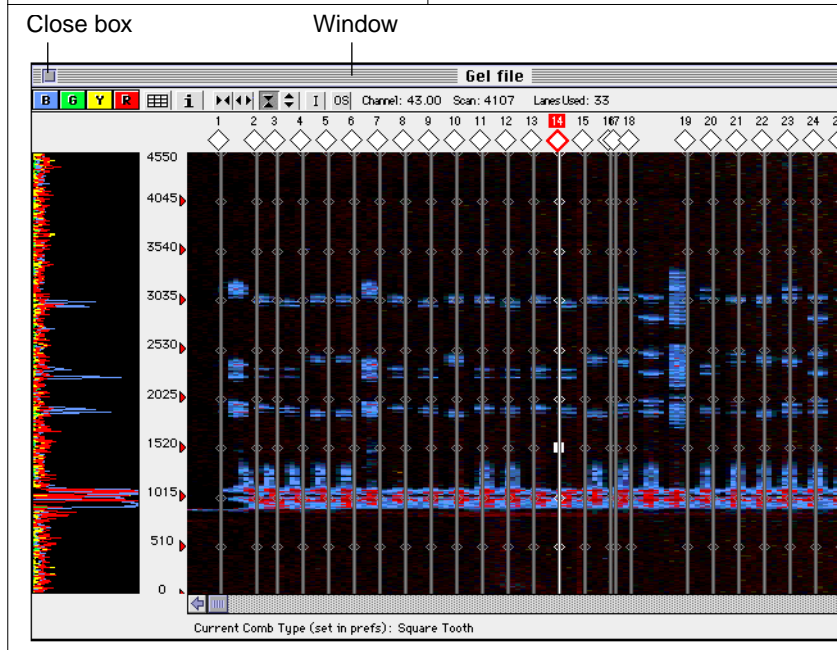
Macintosh Computer Terms Used in This Manual

The following Macintosh® computer terms are used frequently in this manual to describe how to use GeneScan Analysis Software.

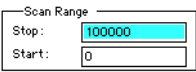
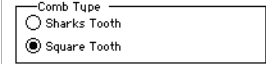

Macintosh computer terms used in this manual.

Item	Description
<p>Dialog Boxes</p> 	<p>Appear when you must make a decision or enter information.</p> <p>All other actions on the monitor screen are suspended until you close the dialog box by clicking a button, such as Cancel, OK, or Done.</p>
<p>Menus</p>	<p>Provide access to various functions you can perform with the software.</p> <p>A heavy arrow (▶) after a menu item indicates that a submenu appears when you click that choice and hold down the mouse button.</p> <p>When you choose a menu command followed by an ellipsis (...) a dialog box appears.</p>
	<p>Menu</p> <p>Submenu</p>
<p>Pop-up menus</p> 	<p>Display a heavy arrow (▶) and are found in dialog boxes.</p> <p>When you click a pop-up menu and hold down the mouse button, a submenu appears.</p> <p>These menus allow you to choose dialog box entries from specific lists of items.</p>

Item	Description
Windows	<p>Display information, and, in some cases, allow you to edit or enter additional information.</p> <p>The top border of an active window always has six horizontal lines and usually has a Close box in the upper-left corner.</p> <p>If many windows are open, click one window to make it active.</p> <p>When a window is active, you can click the top border, hold down the mouse button, and drag the window to another location on the screen.</p> <p>When you are finished working with a window, click the Close box to remove the window from the screen, or click another window.</p>



Macintosh computer terms used in this manual. *(continued)*

Item	Description
<p>Entry fields</p> <p>Image Generation Defaults</p> 	<p>Rectangular areas in which you can enter information.</p> <p>Click in an entry field to display a cursor, and use the keyboard to enter the information.</p>
<p>Checkboxes</p> <p>Auto-Launch Processing</p> <p><input checked="" type="checkbox"/> Auto-Track Gel</p> <p><input checked="" type="checkbox"/> Extract Lanes after Auto-Tracking</p>	<p>Boxes that you click to select certain options in a dialog box.</p> <p>When you click an empty checkbox, an “x” appears in it, indicating that you have selected that option.</p> <p>You can usually select multiple checkboxes.</p>
<p>Radio buttons</p> 	<p>Small circles that appear in front of choices.</p> <p>When you click a radio button with the cursor, a black dot appears in the center of the circle to indicate your choice. You can only select one at a time.</p>
<p>Buttons</p> 	<p>Rectangles with rounded corners that allow you to accept or cancel the contents of a dialog box or perform functions (such as printing) within the dialog box.</p> <p>A button with a heavy outline is the default button that applies if you press the Return key.</p>

Registering the Software

License and Warranty Before you begin, please read Appendix G, "License and Warranty." This appendix explains your rights and responsibilities regarding the software.

Registering Your Software To register your copy of the GeneScan Analysis Software, complete the registration card (included in this software package) and return it to PE Applied Biosystems. Registering the software enables us to send you notification of software updates and any other future information that may be specific to GeneScan Analysis Software owners.

IMPORTANT Your product registration number is located on the Registration card. Be sure to record this number here before you return the Registration card.

Registration Number:

Installing the Software

System Specifications

The following table lists the system specifications¹.

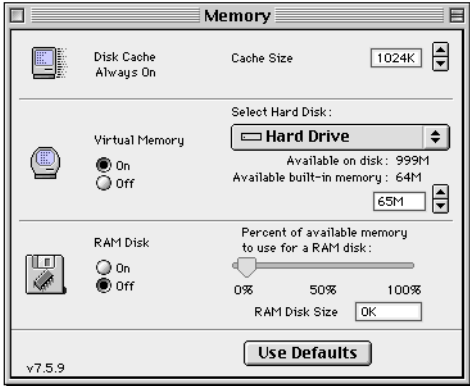
Item	Specifications
CPU	A Power Macintosh® computer (PowerPC CPU). (You will benefit from using the fastest computer available.)
CD-ROM Drive	Any
Operating System	Mac OS version 8.0 with Open Transport 1.1 or later.
Disk Space	A minimum of 15 MB free disk space. Storage requirements depend primarily on the quantity of data to be generated and stored. Sample files are approximately 150—250 KB each and gel files are 20—70 MB each. It is common to store many Sample files on the analysis computer. Gel files are usually stored on the computer that is connected to the instrument and are removed or archived frequently.
Memory (RAM)	The minimum memory requirement is 32 MB total with at least 20 MB of this available to run the GeneScan Analysis Software.
Virtual Memory	Virtual memory MUST be turned on and set to 16 MB. Set the Memory control panel so that the memory available after restart is at least 48M. For information on how, see “Turning On Virtual Memory” on page 1-13. If your computer has 48 MB of physical RAM, virtual memory does not need to be turned on.
Monitor	A 17-inch monitor or larger is recommended. Although a monitor of 640 x 480 resolution can be used, you will benefit from having a monitor of higher resolution.
disk drive	Hard disk with a minimum of 250 MB storage.

1. Call technical support for the latest system specifications.

Item	Specifications
Printer	The following printers are supported: <ul style="list-style-type: none"> ◆ HP DeskJet 1200C/PS ◆ HP DeskJet 1600C/PS
System software	Macintosh computer system software version 8.0.

Turning On Virtual Memory

To turn on Virtual Memory:

Step	Action
1	<p>From the Apple menu (), choose Control Panels and from the list choose Memory. The memory dialog box appears.</p> 
2	<p>Ensure that Virtual Memory is On and that the memory available is at least 48M.</p>

RAM Required to Run the Software

The system requires approximately 32 MB of RAM available to have both the Data Collection software and the analysis programs open simultaneously. This does not include the RAM required by the Macintosh computer operating system.

The RAM required by each program is as follows:

Program	Approximate RAM required
Data Collection software	5.5 MB
GeneScan Analysis Software	32 MB

How to Check the RAM Available

Choose About this Macintosh® from the Apple () menu in the Finder.

Installing the Software

The following procedure describes how to install the GeneScan Analysis Software. Before beginning the installation, backup the copies of the older versions of the GeneScan Analysis Software.

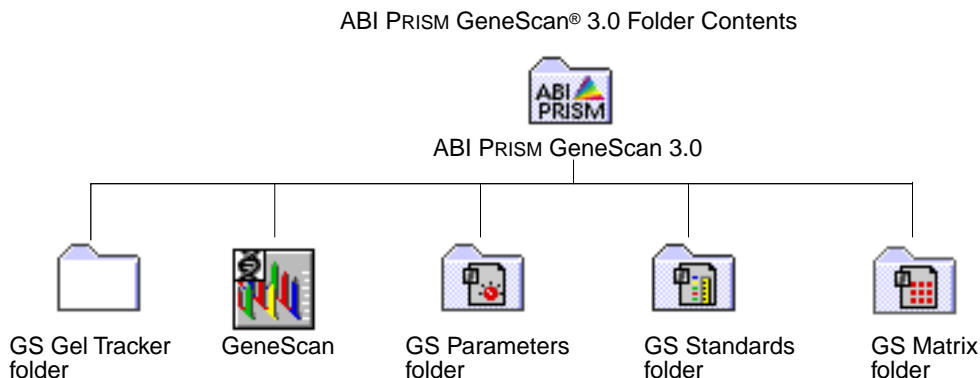
IMPORTANT After installation, be sure to register your software.

To install the GeneScan Analysis Software:

Step	Action
3	Quit all currently running applications.
4	Hold down the Shift key and choose Restart from the Special menu to turn off all extensions.
5	Continue to hold down the shift key until the message "Extension Off" appears.
6	Insert the GeneScan Analysis Software install disk.
7	Double-click the Installer icon. The GeneScan Analysis Software installer appears.
8	Choose a folder in which to install the software.
9	Choose either 310 GeneScan Analysis 3.0 or 373/377/XL GeneScan Analysis 3.0, depending on your instrument.
10	Click Install. Insert other disks as the Installer asks for them.
11	Click Quit at the end of the installation.

Program Files Installed Diagram

The following diagram shows the program files installed on the hard disk.



Contents of Folders

The installation program creates an ABI PRISM GeneScan 3.0 folder on the hard disk containing the following.

Folder	Contents
GS Parameters folder	Use to store user-defined analysis parameters files if you create your own files.
GS Standards folder	Use for storing size standard files when you define them.
GS GelTracker folder	GS Gel Tracker program, tracker settings file, and extensions. You may also see a set of 20 to 30 matlab text files. You can ignore these files; if you throw them away new ones are created when you next track a gel. Note Advise keeping the files because they are useful for troubleshooting auto-tracking problems. Note Do not change the name of the folder, the location of the folder, or the files in the folder.
GS Matrix folder	Use for storing the matrix file generated using the GeneScan Analysis Software. For more information about matrix files, see Chapter 6, "Making a Matrix File."

GeneScan Analysis Software Files

Table of Files The following table lists the files that the GeneScan Analysis Software reads, writes, and, in most cases, creates. The software does not create gel files, and creates Sample files only through lane extraction from a gel. For information on the files installed by the GeneScan Analysis Software, see “Program Files Installed Diagram” on page 1-15.

GeneScan Analysis Software files

File type	Location on		Description
	373XL, ABI PRISM 377, 377XL, 377 96-Lanes, or ABI PRISM 310	ABI 373	
GeneScan Analysis Software	GeneScan 3.0 folder	GeneScan 3.0 folder	The GeneScan Analysis Software analyzes data sent from the ABI PRISM instrument after a run.
The gel file (373 or 377 only)	Run folder created by ABI PRISM 377 Data-Collection during run	Run folder created by ABI 373 Data-Collection during run	<p>A large file created by the Data Collection software.</p> <p>The gel file contains the raw data collected during the instrument run.</p> <p>It also contains the Sample Sheet and run information.</p> <p>For a typical run, this file can be very large. More than 20 MB for 194 channels, more than 40 MB for 388 channels, and more than 50 MB for 96 lanes (up to 70 MB for long runs).</p>

GeneScan Analysis Software files (continued)

File type	Location on		Description
	373XL, ABI PRISM 377, 377XL, 377 96-Lanes, or ABI PRISM 310	ABI 373	
The Project File	Run folder created by the Data Collection software during run	Sample file folder created by GeneScan during run	Organizes, displays, and analyzes Sample files and results. Projects contain references to Sample files. Project files can contain references to Sample files from more than one run.
Sample Files	Run folder created by data-collection software during run	Sample file folder created by GeneScan during run	Created by ABI PRISM 310. For ABI 373 and ABI PRISM 377, GeneScan extracts sample information from the gel file to create Sample files. Each file contains data for a single lane or injection.
Size Standard Files	GS Standards folder or other specified folder	GS Standards folder or other specified folder	Identify peak sizes for specified size standards run under certain conditions. Define the standards after running them on the instrument.
Tracker files and extensions	GS GelTracker folder in the GS 3.0 folder (ABI PRISM 377 and ABI 373 XL)	GS Tracker folder in the GS 3.0 folder	

GeneScan Analysis Software files *(continued)*

File type	Location on		Description
	373XL, ABI PRISM 377, 377XL, 377 96-Lanes, or ABI PRISM 310	ABI 373	
Analysis Parameters	GS Parameters folder or other specified folder	GS Parameters folder or other specified folder	Specify certain ranges and methods used during GeneScan analysis. You can also create and save custom parameters files for future use.
Matrix files	ABI folder, GS Matrix folder or other specified folder	ABI folder, GS Matrix folder or other specified folder	Contains mathematical matrices that correct for spectral overlap. They are applied to data based on the chemistry run.
Analysis Log	GeneScan folder	GeneScan folder	Contains a running record of analysis performed by the GeneScan Analysis Software.
Sample Sheet	GeneScan or specified folder	GeneScan or specified folder	Sample information in spreadsheet form. Used for attaching to gel file.

**Sample File
Database Support**

GeneScan Analysis Software Sample files can be written to and read from the BioLIMS database. The BioLIMS Genetic Information Management System (BioLIMS database) provides a relational database for storage and retrieval of DNA fragment data.

If you have obtained the GeneScan Analysis Software as part of the BioLIMS Client package, you can use the program in either of two modes:

Mode	Description
Sample File mode	<p>Fragment data extracted from gel files is written to individual sample files.</p> <p>Extracted fragment data viewed and processed within the GeneScan Analysis Software is read from and saved to sample files.</p>
BioLIMS mode	<p>Fragment data extracted from gel files is written directly to a BioLIMS database that resides on a UNIX Server.</p> <p>Extracted fragment data viewed and processed within the GeneScan Analysis Software is read from and written back to the same BioLIMS database.</p> <p>For more information on using the BioLIMS database, see Appendix E, "Using the BioLIMS Database."</p>

Technical Support

To Reach Us On the Web

Our web site address is <http://www.perkin-elmer.com/ab>

We strongly encourage you to visit our web site for answers to frequently asked questions, and to learn more about our products. You can also order technical documents and/or an index of available documents and have them faxed to you through our site (see the "Fax-on-Demand" section below).

Hours for Telephone Technical Support

In the United States and Canada, technical support is available between 5:30 a.m. and 5:00 p.m. Pacific Time.

These hours are for all products except the following:

Product	Hours
Chemiluminescence	9:00 a.m. to 5:00 p.m. Eastern Time
LC/MS	9:00 a.m. to 5:00 p.m. Pacific Time

See the "Regional Offices Sales and Service" section below for how to contact local service representatives outside of the United States and Canada.

To Reach Us by Telephone or Fax in North America

Call Technical Support at 1-800-831-6844, and select the appropriate option (below) for support on the product of your choice at any time during the call. (To open a service call for other support needs, or in case of an emergency, press 1 after dialing 1-800-831-6844.)

For Support On This Product	Dial 1-800-831-6844, and...	
DNA Synthesis	Press	FAX
	21	650-638-5981
Fluorescent DNA Sequencing	Press	FAX
	22	650-638-5891
Fluorescent Fragment Analysis (includes GeneScan applications)	Press	FAX
	23	650-638-5891

For Support On This Product	Dial 1-800-831-6844, and...					
BioLIMS	<table border="1"> <thead> <tr> <th data-bbox="749 172 995 220">Press</th> <th data-bbox="995 172 1243 220">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 220 995 269">25</td> <td data-bbox="995 220 1243 269">650-638-5891</td> </tr> </tbody> </table>	Press	FAX	25	650-638-5891	
Press	FAX					
25	650-638-5891					
Integrated Thermal Cyclers	<table border="1"> <thead> <tr> <th data-bbox="749 285 995 334">Press</th> <th data-bbox="995 285 1243 334">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 334 995 383">24</td> <td data-bbox="995 334 1243 383">650-638-5891</td> </tr> </tbody> </table>	Press	FAX	24	650-638-5891	
Press	FAX					
24	650-638-5891					
PCR and Sequence Detection	<table border="1"> <thead> <tr> <th data-bbox="749 399 995 448">Press</th> <th data-bbox="995 399 1243 448">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 448 995 651">5, or call 1-800-762-4001, and press 1 for PCR, or 2 for Sequence Detection</td> <td data-bbox="995 448 1243 651">203-761-2542</td> </tr> </tbody> </table>	Press	FAX	5, or call 1-800-762-4001, and press 1 for PCR, or 2 for Sequence Detection	203-761-2542	
Press	FAX					
5, or call 1-800-762-4001, and press 1 for PCR, or 2 for Sequence Detection	203-761-2542					
Peptide Synthesis	<table border="1"> <thead> <tr> <th data-bbox="749 667 995 716">Press</th> <th data-bbox="995 667 1243 716">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 716 995 764">31</td> <td data-bbox="995 716 1243 764">650-638-5981</td> </tr> </tbody> </table>	Press	FAX	31	650-638-5981	
Press	FAX					
31	650-638-5981					
Protein Sequencing	<table border="1"> <thead> <tr> <th data-bbox="749 781 995 829">Press</th> <th data-bbox="995 781 1243 829">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 829 995 878">32</td> <td data-bbox="995 829 1243 878">650-638-5981</td> </tr> </tbody> </table>	Press	FAX	32	650-638-5981	
Press	FAX					
32	650-638-5981					
Chemiluminescence	<table border="1"> <thead> <tr> <th data-bbox="749 894 995 943">Telephone</th> <th data-bbox="995 894 1243 943">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 943 995 1105">1-800-542-2369 (U.S. only), or 1-617-271-0045 (Tropix)</td> <td data-bbox="995 943 1243 1105">617-275-8581 (Tropix) 9:00 a.m. to 5:00 p.m. ET</td> </tr> </tbody> </table>	Telephone	FAX	1-800-542-2369 (U.S. only), or 1-617-271-0045 (Tropix)	617-275-8581 (Tropix) 9:00 a.m. to 5:00 p.m. ET	
Telephone	FAX					
1-800-542-2369 (U.S. only), or 1-617-271-0045 (Tropix)	617-275-8581 (Tropix) 9:00 a.m. to 5:00 p.m. ET					
LC/MS	<table border="1"> <thead> <tr> <th data-bbox="749 1122 995 1170">Telephone</th> <th data-bbox="995 1122 1243 1170">FAX</th> </tr> </thead> <tbody> <tr> <td data-bbox="749 1170 995 1294">1-800-952-4716</td> <td data-bbox="995 1170 1243 1294">650-638-6223 9:00 a.m. to 5:00 p.m. PT</td> </tr> </tbody> </table>	Telephone	FAX	1-800-952-4716	650-638-6223 9:00 a.m. to 5:00 p.m. PT	
Telephone	FAX					
1-800-952-4716	650-638-6223 9:00 a.m. to 5:00 p.m. PT					

Fax-on-Demand Free 24-hour access to PE Applied Biosystems technical documents is available by fax.

You can access Fax-on-Demand documents through the internet or by telephone:

If you want to order...	Then...
through the internet	Use http://www.perkin-elmer.com/ab/techsupp You can search for documents to order using keywords. Up to five documents can be faxed to you if you already know the titles.
by phone from the United States or Canada	<ol style="list-style-type: none">Call 1-800-487-6809 from a touch-tone phone. Have your fax number ready.Press 1 to order an index of available documents and have it faxed to you. Each document in the index has an ID number. (Use this as your order number in step "d" below.)Call 1-800-487-6809 from a touch-tone phone a second time.Press 2 to order up to five documents and have them faxed to you.
by phone from outside the United States or Canada	<ol style="list-style-type: none">Dial your international access code, then 1-650-596-4419, from a touch-tone phone. Have your complete fax number and country code ready (011 precedes the country code).Press 1 to order an index of available documents and have it faxed to you. Each document in the index has an ID number. (Use this as your order number in step "d" below.)Call 1-650-596-4419 from a touch-tone phone a second time.Press 2 to order up to five documents and have them faxed to you.

**To Reach Us by
E-Mail**

Contact technical support by e-mail for help in the following product areas.

For this product area	Use this e-mail address
Chemiluminescence	info@tropix.com
Genetic Analysis	galab@perkin-elmer.com
LC/MS	apisupport@sciex.com
PCR and Sequence Detection	pcriab@perkin-elmer.com
Protein Sequencing, Peptide and DNA Synthesis	corelab@perkin-elmer.com

**Regional Offices
Sales and Service**

If you are outside the United States and Canada, you should contact your local PE Applied Biosystems service representative.

The Americas	
United States PE Applied Biosystems 850 Lincoln Centre Drive Foster City, California 94404 Tel: (650) 570-6667 (800) 345-5224 Fax: (650) 572-2743	Canada (Mississauga, Ontario) Tel: (905) 821-8183 (800) 668-6913 Fax: (905) 821-8246 Latin America (Del.A. Obregon, Mexico) Tel: (305) 670-5310 Fax: (305) 670-4349

Europe	
Austria (Wien) Tel: 01 602 3101 Fax: 01 602 5174	Hungary (Budapest) Tel: 1 251 11 16 Fax: 1 251 14 61
Benelux (Nieuwerkerk a/d IJssel, Netherlands) Tel: 31(0) 180 331400 Fax: 31(0) 180 331409	Italy (Milano) Tel: (039) 23831 Fax: (039) 2383490
Chekia Rep. (Praha) Tel: 2 61 22 21 64 Fax: 2 61 22 21 68	Norway (Oslo) Tel: (0) 22 02 1500 Fax: (0) 22 02 1501

Europe <i>(continued)</i>	
Denmark (Allerød) Tel: 48 100 400 Fax: 48 100 401	Poland (Warszawa) Tel: (48 22) 866 40 10 Fax: (48 22) 866 40 20
Finland (Espoo) Tel: 09 751 72 700 Fax: 09 751 72 701	Russia (Moskva) Tel: 095 935 8888 Fax: 095 564 8787
France (Paris) Tel: (1) 69 59 85 85 Fax: (1) 69 59 85 00	South Africa (Johannesburg) Tel: 27 11 478 0411 Fax: 27 11 478 0349
Germany (Weiterstadt) Tel: (0) 6150/ 101 0 Fax: (0) 6150/ 101 101	Spain (Madrid) Tel: (1) 806 1200 Fax: (1) 804 0414
Sweden (Sundbyberg) Tel: (0)8 619 4400 Fax: (0)8 619 4401	United Kingdom (Warrington, Cheshire) Tel: (01925) 825650 Fax: (01925) 282502
Switzerland (Rotkreuz) Tel: (0) 41 799 7708 Fax: (0) 41 790 0676	All Other European Countries, Middle East/West Asia, Africa (Langen, Germany) Tel: 49 6103 708 301 Fax: 49 6103 708 310

Japan
Japan (Chiba) Tel: (0473) 80 8500 Fax: (0473) 80 8505

Eastern Asia, China, Oceania	
Australia (Scoresby, Victoria) Tel: (03) 9212 8585 Fax: (03) 9212 8502	Malaysia (Kuala Lumpur) Tel: 60 3 758 1118 Fax: 60 3 754 9043
China (Beijing) Tel: 86 10 6238 1156 Fax: 86 10 6238 1162	Singapore Tel: 65 336 0322 Fax: 65 338 3991

Eastern Asia, China, Oceania	
Hong Kong Tel: 852 2756 6928 Fax: 852 2756 6968	Taiwan (Taipei Hsish) Tel: 886 2 698 3505 Fax: 886 2 698 3405
Korea (Seoul) Tel: 822 592 7238 Fax: 822 532 4908	Thailand (Bangkok) Tel: 662 719 6406 Fax: 662 319 9788

How to Process and Edit the Gel File

2

Introduction

In This Chapter This chapter contains information about how to view and edit the gel file, and how to generate Sample files or BioLIMS™ database records after editing the gel file.

Note This section does not apply if you are analyzing Sample files generated on the ABI PRISM® 310 instrument.

IMPORTANT Refer to Appendix A, “Using GeneScan with the ABI 373,” before using the GeneScan program to analyze data collected on the ABI 373.

Topics in This Chapter Topics in this chapter include the following:

Topic	See page
About Gel Files	2-2
How to Set Gel Processing Preferences	2-3
How to Display the Gel File	2-11
About the Gel File Window	2-12
How to Adjust the Gel Image	2-18
Verifying Gel Information	2-26
Displaying the Sample Sheet	2-29
About the Sample Sheet	2-31
Working with the Sample Sheet	2-34
About Displaying Regions of Off-Scale Data	2-38
Working with Lane Markers	2-43
Marking and Unmarking Lanes for Extraction	2-50
Working with Tracker Lines	2-53
Interpolating Tracker Lines	2-57
Tracking Lanes and Extracting Data	2-59

Topic	See page
How the GeneScan Analysis Software Names Sample Files	2-67
Saving Gel Files After Editing Tracking	2-68

About Gel Files

Instruments Producing Gel File

The gel file, which is generated only by the ABI™ 373 DNA Sequencer instrument, the ABI 373 with XL upgrade, the ABI PRISM® 377 DNA Sequencer instrument, and the ABI PRISM® 377 with XL upgrade, stores the raw data collected during the entire run of the instrument. A gel file can be 10 MB to 60 MB.

96-Lane Gels Capability

The GeneScan® Analysis Software can open and analyze 96-lane gels.

Gel File Contents

Initially, the file contains the following:

- ◆ Raw data collected during the run.
- ◆ Gel image, which is similar to an autoradiogram, but in color.
- ◆ Copy of the Data Collection software sample sheet.
- ◆ Copy of the matrix file.

After lane tracking and editing, the file also contains the lane tracking information and any changes you make to the original information in the file.

How GeneScan Tracks a Gel File

When the GeneScan Analysis Software tracks the gel file, it creates a tracker line for each lane in the gel. Each tracker line is trying to find the center of the bands, not the brightest part of the band. The tracker will only track if a matrix is attached to the gel file.

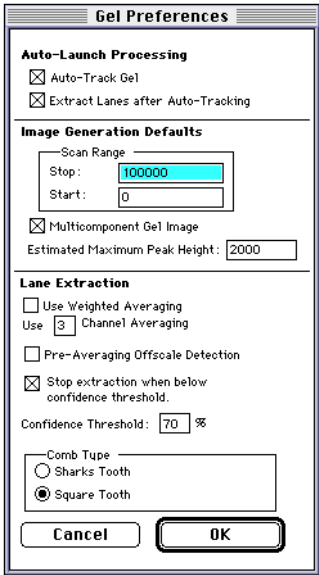
During data extraction, the software generates a Sample file for each tracked lane by averaging the data from the tracked channel and the number of channels you specify on either side of it. The software also copies to each new Sample file all the information required to identify and analyze the sample.

How to Set Gel Processing Preferences

Introduction Use the Gel Processing dialog box to define the parameter values to use for gel file processing. The following procedure applies to the ABI 373 and the ABI PRISM 377 instruments.

Displaying the Gel Preferences Dialog Box

To display the Gel Preferences dialog box:

Step	Action
1	If the GeneScan Analysis Software is not running, then double-click the program icon.
2	<p>Choose Gel Preferences from the Settings menu.</p> <p>The Gel Preferences dialog box appears.</p>  <p>There are three Gel Preferences parameters:</p> <ul style="list-style-type: none"> ◆ Auto-Launch Processing. ◆ Image Generation Defaults. ◆ Lane Extraction.

Auto-Launch Processing

Select or de-select the checkboxes to specify whether or not the gel is to be automatically tracked and extracted at the end of a data collection run.

IMPORTANT Before auto-tracking the first time, you need to verify that the comb type is set correctly. The default comb type is Square.

Note Select both checkboxes to analyze samples automatically after data-collection.

Image Generation Defaults

In the Image Generation Defaults options enter values that determine how the GeneScan Analysis Software processes the gel file when it generates the initial gel image from data collected after electrophoresis.

Image Generation options:

Item	Description						
Scan Range	<p>Enter the scan numbers at which you want to display the gel image. Use a trial run to determine the scan range since the time frame in which fragments are detected may vary from run to run.</p> <p>During analysis, the GeneScan Analysis Software ignores all scans outside the specified scan range, and only includes the scans with the designated ranges when it generates Sample files.</p> <p>See "Regenerating the Gel Image" on page 2-23 to change the number of scans displayed.</p>						
	<p>Determining the Start and Stop Scan Numbers:</p> <p>a. Determine the start number by multiplying the applicable number from the table below by the number of hours it takes for the peaks to appear.</p> <table border="1"><thead><tr><th>Instrument & Type of Run</th><th>Scans per hour</th></tr></thead><tbody><tr><td>ABI 373</td><td>600</td></tr><tr><td>ABI PRISM 377 and ABI™ 373 with XL upgrade</td><td>1,200 to 2,400^a</td></tr></tbody></table> <p>a. Depending on which module is used during data collection</p> <p>b. Determine the stop number by multiplying the applicable number from the above table by the number of hours that pass before the last peak is detected.</p>	Instrument & Type of Run	Scans per hour	ABI 373	600	ABI PRISM 377 and ABI™ 373 with XL upgrade	1,200 to 2,400 ^a
Instrument & Type of Run	Scans per hour						
ABI 373	600						
ABI PRISM 377 and ABI™ 373 with XL upgrade	1,200 to 2,400 ^a						

Image Generation options: *(continued)*

Item	Description
Multicomponent Gel Image	<p>Note This option affects only the gel image.</p> <p>Select this checkbox to use the spectral separation algorithm in conjunction with a predefined matrix file to resolve the spectral overlap between dyes.</p> <p>Matrix files are specific to a particular set of run conditions and to the instrument on which they are created.</p> <p>In other words, you must create a different matrix file for each dye set, gel type, and set of running conditions you use.</p> <p>For more information, see Chapter 6, “Making a Matrix File.”</p> <p>IMPORTANT If the gel does not have an associated matrix, you must attach a matrix to the gel, or the software will not multicomponent the gel image, even if you select this option. The tracker relies upon multicomponent images for accurate tracking.</p> <p>IMPORTANT To attach a matrix, see “Installing New Matrix Information” on page 2-24.</p> <p>Note You must always attach a matrix to the ABI 373 Collection file in order to multicomponent the gel image. ABI 672 Data Collection software is not capable of pre-installing a matrix.</p> <p>See Appendix A, “Using GeneScan with the ABI 373.”</p>

Image Generation options: *(continued)*

Item	Description
<p>Estimated Maximum Peak Height</p>	<p>In this text box, enter the maximum signal level you expect from samples in the run. This can be an approximate number, based on your typical run conditions and samples.</p> <p>The default is 2000.</p> <p>Note This option affects only the appearance of the gel image (not the raw data) when the image is generated the first time the gel file is opened. Use the Regenerate Gel Image command to change the appearance of an open gel window (See page 2-23).</p> <p>The gel image can also be modified using the Adjust Gel Contrast command; see “Adjusting the Contrast” on page 2-20.</p> <p>The Estimated Maximum Peak Height value:</p> <ul style="list-style-type: none"> ◆ Affects the color brightness of the DNA fragment bands. <p>All bands with a data value at or above the Estimated Maximum Peak Height value are assigned the brightest dye color.</p> <p>The dye colors for bands with values below that level are dimmed proportionally, unless multicomponented.</p> <p>In general, the lower you set this value, the brighter the bands appear in the Gel File window. A value of 1000 is satisfactory for most gel files. If the gel image is very dim, try 500; if it is too bright, try 2000.</p> ◆ Determines the scale of the peaks in the Slice view of the Gel File window. <p>For the highest quality gel image, the highest sample peaks (<i>not</i> the Primer peak) should just reach the top of the scale in the Slice view.</p> <p>If you find that many of the peaks are cut off, you may want to readjust the Estimated Peak Height value to a higher number.</p>

Lane Extraction The following table lists the options in the Lane Extraction section.

Item	Description
Use Weighted Averaging checkbox	<p>Weighted channel averaging is a new feature in GeneScan v. 3.0.</p> <p>Weighted averaging is now possible because the new Tracker interface allows tracker line placement to within a tenth of a channel.</p>

No Weighted Averaging

If the Use Weighted Averaging box is not checked, data averaging is done per channel the same as with earlier versions of the GeneScan Analysis Software.

For example, if the tracker line falls within channel 10, and three-channel averaging is set:

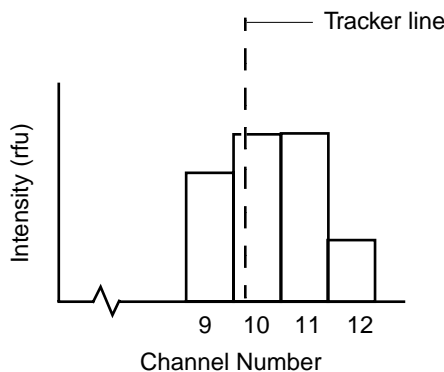
$$\text{channel average} = \frac{I_{\text{ch. 9}} + I_{\text{ch. 10}} + I_{\text{ch. 11}}}{3}$$

Where I is the intensity for a given channel and scan number.

Weighted Averaging

If the Use Weighted Averaging box is selected, the tracker line falls 20% into channel 10 (see diagram below), and two-channel averaging is set:

$$\text{channel average} = \frac{(I_{\text{ch. 9}} \times 0.7) + (I_{\text{ch. 10}} \times 1) + (I_{\text{ch. 11}} \times 1) + (I_{\text{ch. 12}} \times 0.3)}{3}$$



Item	Description
Use Channel Averaging text entry field	<p data-bbox="615 142 1180 256">Averaging reduces the amount of noise in the sample file and allows for more accurate representation of the band as a peak. (See “Weighted Averaging” on page 2-7).</p> <p data-bbox="615 272 1180 354">The number of channels to be averaged for each lane when extracting data from the gel file is normally set to 3.</p> <p data-bbox="615 378 1180 459">Each tracker line in the Gel File window marks the channel where the GeneScan Analysis Software located the strongest fluorescent signal for that lane.</p> <p data-bbox="615 483 1180 597">If you use the default three-channel average (without weighted-averaging), the raw data in each sample file is an average of the data in the channel marked by the tracker line and one channel on either side of it.</p> <p data-bbox="615 621 1180 813">IMPORTANT When using multiple-channel averaging, ensure the tracker lines mark central channels of the bands. If they mark channels at the right or left edge of the bands, the empty channel between lanes is included in the average or the signal from an adjacent lane could be included, causing an erroneous value.</p>

Item	Description						
	<p data-bbox="662 139 1026 167">Altering the Channel Averaging</p> <table border="1" data-bbox="662 207 1239 898"> <thead> <tr> <th data-bbox="662 207 951 245">If you choose...</th> <th data-bbox="951 207 1239 245">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="662 245 951 594">two-channel averaging</td> <td data-bbox="951 245 1239 594"> <p data-bbox="964 253 1225 451">data is taken from the tracked channel and the channel to the right of it (without weighted averaging). You can include data from up to nine channels.</p> <p data-bbox="964 475 1214 586">Three-channel averaging is recommended for most applications.</p> </td> </tr> <tr> <td data-bbox="662 594 951 898">one-channel averaging (no averaging)</td> <td data-bbox="951 594 1239 898"> <p data-bbox="964 602 1220 886">if the gel bands are severely tilted. For example, if the left lane of the gel ran faster than the right, a better result would be obtained by taking the center channel alone, rather than averaging three channels.</p> </td> </tr> </tbody> </table>	If you choose...	Then...	two-channel averaging	<p data-bbox="964 253 1225 451">data is taken from the tracked channel and the channel to the right of it (without weighted averaging). You can include data from up to nine channels.</p> <p data-bbox="964 475 1214 586">Three-channel averaging is recommended for most applications.</p>	one-channel averaging (no averaging)	<p data-bbox="964 602 1220 886">if the gel bands are severely tilted. For example, if the left lane of the gel ran faster than the right, a better result would be obtained by taking the center channel alone, rather than averaging three channels.</p>
If you choose...	Then...						
two-channel averaging	<p data-bbox="964 253 1225 451">data is taken from the tracked channel and the channel to the right of it (without weighted averaging). You can include data from up to nine channels.</p> <p data-bbox="964 475 1214 586">Three-channel averaging is recommended for most applications.</p>						
one-channel averaging (no averaging)	<p data-bbox="964 602 1220 886">if the gel bands are severely tilted. For example, if the left lane of the gel ran faster than the right, a better result would be obtained by taking the center channel alone, rather than averaging three channels.</p>						
Pre-Averaging Offscale Detection checkbox	<p data-bbox="662 914 1231 1052">When you select this checkbox, if any of the channels that are used for averaging has a value of 8191 relative fluorescent units (rfu), then no averaging takes place and the value of 8191 rfu is written for that scan of that lane.</p> <p data-bbox="662 1076 1228 1187">Note If you extracted a sample file with the Pre-Averaging Offscale Detection checkbox selected, then pre-averaging has a value of zero in the Sample Info View.</p> <p data-bbox="662 1211 1224 1268">For information about off-scale data, see “About Displaying Regions of Off-Scale Data” on page 2-38.</p> <p data-bbox="662 1292 1080 1312">The default is checkbox is not selected.</p>						

Item	Description
Stop extraction when below confidence value checkbox	<p>If the Stop extraction when below confidence threshold box is checked, the lane extraction will not be carried out and a warning dialog box will appear when the lane assignment confidence level is below that specified in the Confidence Threshold.</p> <p>The dialog box gives you the option to cancel or continue the gel file extraction and analysis.</p>
Confidence Threshold text entry field	<p>After a gel is auto-tracked, a lane assignment confidence value is written to the Analysis Log.</p> <p>This value indicates the Tracker's confidence in how well the assigned lanes match the sample sheet. (This value is not an indication of how well the tracker lines follow the fluorescence intensity within the lanes.)</p> <p>Lane assignment confidence values tend to be extreme numbers: very low or very high.</p> <p>Although, a value of 70 or more generally indicates that the lane assignment for the gel is correct, it is recommended that you check the tracker lane assignment anytime the reported lane assignment confidence value is less than 100%.</p>

Comb Type **IMPORTANT** Before auto-tracking the first time, you need to verify that the comb type is set correctly. The default comb type is Square.

The Neural Net Tracker uses special tracker setting files that are optimized according to the number of channels and lanes in the gel file and the comb type.

It is important that you set the correct comb type for the gel so that the Tracker applies the correct tracker setting files.

This value appears at the bottom of the gel window (see "Gel File Window Diagram" on page 2-13).

IMPORTANT The proper comb type needs to be selected in order to ensure proper tracker performance.

How to Display the Gel File

Introduction Use the gel image to observe sample migration, lane tracking, and signal strength in the gel image. You can also use the window to adjust the image to improve visibility, re-align individual lane markers, or edit the position of the tracker lines.

Displaying the Gel File You can display a gel file as follows:

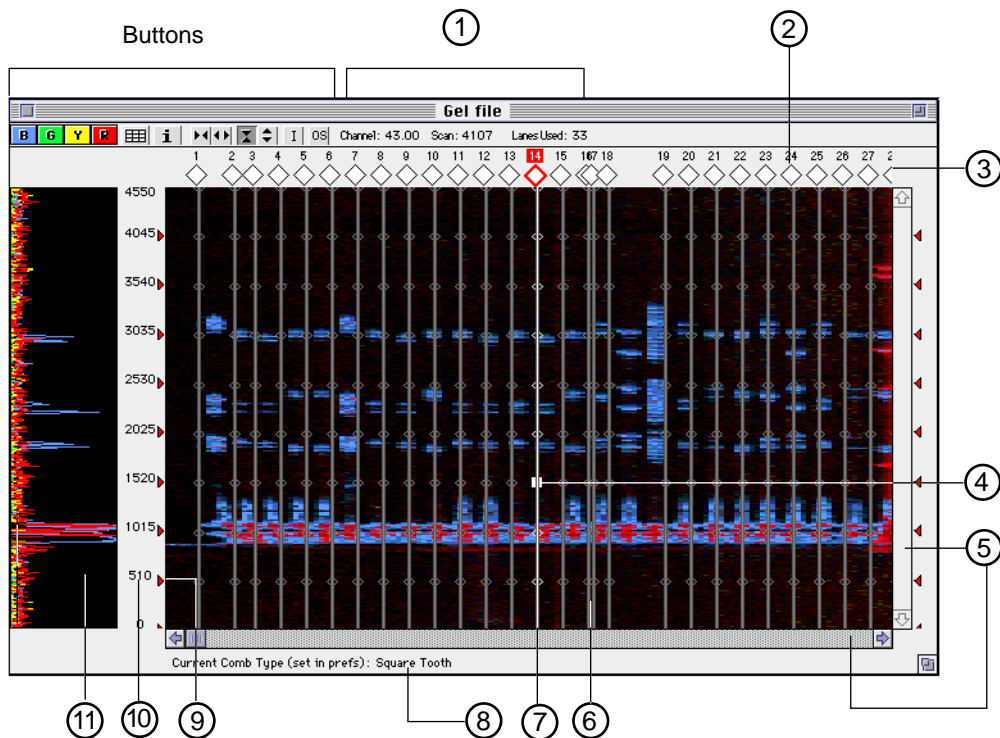
To display a gel file...	Then...
automatically	after automatic analysis, the Gel File window opens and displays the newly created gel file. For more information, see "Setting Up for Automatic Analysis" on page 4-4.
manually	you can either: Double-click the Gel file, or a. Choose Open from the File menu. The Open Existing dialog box appears. b. Click the Collection Gel icon. A directory dialog box appears. c. Locate and select the desired gel file and choose Open.

About the Gel File Window

Differences Between Gel File Windows The following table explains the differences between the GeneScan Analysis Software Gel File window and the Gel File window displayed by the Data Collection software during a run.








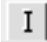
Application	Shows	New Data	Gel Image
Data Collection software Gel window	Real-time data as it is being collected.	New data appears at the bottom of the screen as it is collected, so the top of the screen shows the start of the run.	Gel image is not baselined or multicomponented
GeneScan Gel File window	An image of the gel after the Data Collection software is finished.	This image is inverted. The bottom of the window displays the start of the run, that is, the smallest fragments appear at the bottom of the window, just as they would on an autoradiogram.	Gel image is baselined and may or may not be multicomponented.

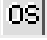
Gel File Window Diagram



Gel File Window Buttons

The following table describes the buttons on the Gel File window.

Window tool	Description
	Color buttons—Turns on/off the display of one or more colors in the gel image and slice view (see also page 2-18).
	<p>Sample Sheet button—Displays a copy of the Data Collection Sample Sheet associated with the gel file.</p> <p>You can use the sample sheet information to check that the lanes are correctly labeled for the gel image.</p> <p>For more information, see “Displaying the Sample Sheet” on page 2-29.</p>
	<p>Gel Info button—Displays the Gel Info dialog box.</p> <p>For more information, see “Gel Info Window” on page 2-26.</p>
	Horizontal Shrink button—Compresses the gel image horizontally in order to see all the gel lanes in a standard-sized window.
	<p>Horizontal Expand button—Expands the gel image horizontally in order to adjust the tracker lanes.</p> <p>The four levels of horizontal zoom are: 1x, 2x, 4x, and 8x.</p>
	Vertical Shrink button—Returns the vertical scale to normal after using the Vertical Expand button.
	<p>Vertical Expand button—Expands the gel image vertically to see the quality of the peaks in Slice View.</p> <p>There are two levels of vertical zoom: full scale and 600 scans.</p>
	<p>Interpolation button—Places the user in tracker-line interpolation mode.</p> <p>For more information, see “Interpolating Tracker Lines” on page 2-57.</p>

Window tool	Description
	<p>Off-Scale button—Displays off-scale data.</p> <p>Turns off all the colors in the gel image and displays in white any data points that are off-scale.</p> <p>Off-Scale data is outside the range of the detector (8191 relative fluorescent units and greater.)</p> <p>Note To see all off-scale data points, the Horizontal Expand and the Vertical Expand buttons must be used.</p> <p>For more information, see “About Displaying Regions of Off-Scale Data” on page 2-38.</p>

Gel File Window Described

The following table describes the diagram on page 2-13.

Gel File window described.

No.	Description						
1	<table border="1"> <tbody> <tr> <td style="text-align: center;">Channel</td> <td> <p>Current cursor position on the horizontal scale and is displayed in sub-channel increments.</p> <p>Channels are the divisions across the read region of a gel where the Data Collection software samples the data.</p> <p>The number of available channels depends on the instrument and run mode used.</p> </td> </tr> <tr> <td style="text-align: center;">Scan</td> <td> <p>Current cursor position on the vertical scale.</p> <p>Note These numbers change as you move the cursor over the image.</p> </td> </tr> <tr> <td style="text-align: center;">Lane Used</td> <td> <p>Number of lanes used in the gel.</p> </td> </tr> </tbody> </table>	Channel	<p>Current cursor position on the horizontal scale and is displayed in sub-channel increments.</p> <p>Channels are the divisions across the read region of a gel where the Data Collection software samples the data.</p> <p>The number of available channels depends on the instrument and run mode used.</p>	Scan	<p>Current cursor position on the vertical scale.</p> <p>Note These numbers change as you move the cursor over the image.</p>	Lane Used	<p>Number of lanes used in the gel.</p>
Channel	<p>Current cursor position on the horizontal scale and is displayed in sub-channel increments.</p> <p>Channels are the divisions across the read region of a gel where the Data Collection software samples the data.</p> <p>The number of available channels depends on the instrument and run mode used.</p>						
Scan	<p>Current cursor position on the vertical scale.</p> <p>Note These numbers change as you move the cursor over the image.</p>						
Lane Used	<p>Number of lanes used in the gel.</p>						
2	<p>Lane numbers—Numbers across the top of the gel image show the lane number currently assigned to each lane on the gel.</p>						

Gel File window described. (continued)

No.	Description												
3	<p>Lane markers—Diamond shaped markers (◆) show the current status of each lane.</p> <p>The following table lists the status values:</p> <table border="1" data-bbox="465 324 1184 784"> <thead> <tr> <th data-bbox="465 324 610 391">Lane color</th> <th data-bbox="610 324 1184 391">Description</th> </tr> </thead> <tbody> <tr> <td data-bbox="465 391 610 430">White</td> <td data-bbox="610 391 1184 430">Lane used and marked for extraction.</td> </tr> <tr> <td data-bbox="465 430 610 469">Blue</td> <td data-bbox="610 430 1184 469">Lane is used but not marked for extraction.</td> </tr> <tr> <td data-bbox="465 469 610 540">Yellow</td> <td data-bbox="610 469 1184 540">Lane was edited and extracted, but the gel file was not saved with new information.</td> </tr> <tr> <td data-bbox="465 540 610 638">Gray</td> <td data-bbox="610 540 1184 638">The tracker does not expect to find a lane here, if it does, the tracker will confuse it and lane assignment confidence will be low.</td> </tr> <tr> <td data-bbox="465 638 610 784">Orange</td> <td data-bbox="610 638 1184 784"> <p>Lane was inferred by the Tracker software.</p> <p>If you move or reshape an inferred lane tracker line, it ceases to be inferred and the orange border is lost.</p> </td> </tr> </tbody> </table>	Lane color	Description	White	Lane used and marked for extraction.	Blue	Lane is used but not marked for extraction.	Yellow	Lane was edited and extracted, but the gel file was not saved with new information.	Gray	The tracker does not expect to find a lane here, if it does, the tracker will confuse it and lane assignment confidence will be low.	Orange	<p>Lane was inferred by the Tracker software.</p> <p>If you move or reshape an inferred lane tracker line, it ceases to be inferred and the orange border is lost.</p>
Lane color	Description												
White	Lane used and marked for extraction.												
Blue	Lane is used but not marked for extraction.												
Yellow	Lane was edited and extracted, but the gel file was not saved with new information.												
Gray	The tracker does not expect to find a lane here, if it does, the tracker will confuse it and lane assignment confidence will be low.												
Orange	<p>Lane was inferred by the Tracker software.</p> <p>If you move or reshape an inferred lane tracker line, it ceases to be inferred and the orange border is lost.</p>												
4	<p>Control points—Use to control shape and position the splines.</p> <p>For more information, see “Using Control Points” on page 2-45.</p>												
5	<p>Scroll bars—Use to scroll the gel view horizontally or vertically.</p>												
6	<p>Gel Image—An image representing a time history of all fluorescence detected during the run.</p> <ul style="list-style-type: none"> ◆ Each peak is shown as a brightly colored band within the sample lane. ◆ Each position on the gel image is defined by a scan number and a channel number. 												
7	<p>Tracker line for the selected lane.</p>												
8	<p>Current Comb Type—Displays the current comb type that is set in the Gel preferences.</p> <p>For more information, see “Comb Type” on page 2-10.</p>												
9	<p>Control row</p> <p>For information on adding and deleting a control row, see “Adding and Deleting Control Rows” on page 2-46.</p>												

Gel File window described. *(continued)*

No.	Description						
10	<p>Scan number—Reflects a particular point in time.</p> <p>The address of a particular data point on the Gel file image.</p>						
	<table border="1"> <thead> <tr> <th data-bbox="517 293 709 332">This run...</th> <th data-bbox="709 293 1245 332">Collects...</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 332 709 373">2400</td> <td data-bbox="709 332 1245 373">approximately 2250 scans per hour</td> </tr> <tr> <td data-bbox="517 373 709 414">1200</td> <td data-bbox="709 373 1245 414">approximately 1125 scans per hour</td> </tr> </tbody> </table>	This run...	Collects...	2400	approximately 2250 scans per hour	1200	approximately 1125 scans per hour
	This run...	Collects...					
	2400	approximately 2250 scans per hour					
1200	approximately 1125 scans per hour						
<table border="1"> <tbody> <tr> <td data-bbox="709 332 1245 373">approximately 2250 scans per hour</td> </tr> </tbody> </table>	approximately 2250 scans per hour						
approximately 2250 scans per hour							
<table border="1"> <tbody> <tr> <td data-bbox="709 373 1245 414">approximately 1125 scans per hour</td> </tr> </tbody> </table>	approximately 1125 scans per hour						
approximately 1125 scans per hour							
11	<p>Slice view—Graphical view of the data values in the tracked channel of the selected lane.</p> <ul style="list-style-type: none"> ◆ The display changes as you move the tracker line from one channel to another in the lane. ◆ Each peak in the Slice view corresponds to a band in the gel image and indicates a DNA fragment. ◆ Although these bands and peaks do not represent analyzed data, they provide an overview of the relative signal intensity between the bands in that lane, and allow you to make a qualitative evaluation of the run. ◆ The Slice view is empty (black) when no lane is selected or when the lane has no data. 						

How to Adjust the Gel Image

Introduction Although the data shown in the gel image is not analyzed, the displayed information allows you to evaluate the quality of the run.

You can adjust the appearance of the gel image as follows.

- ◆ Displaying/Hiding a Dye Color.
- ◆ Changing Dye Indicators.
- ◆ Adjusting the Contrast.
- ◆ Regenerating the Gel Image using a different scan number range, estimated maximum peak height, and multicomponenting option.
- ◆ Installing New Matrix Information.


Note Changing any of the above features in the Gel File window does not change the raw data contained in the gel file.

Displaying/Hiding a Dye Color

Introduction

You can control the display of specific fluorescent dye colors. For example, if you want to display only the bands labeled with blue dye, you can suppress the display of all green, yellow, or red bands.

How to Display or Hide a Dye Color

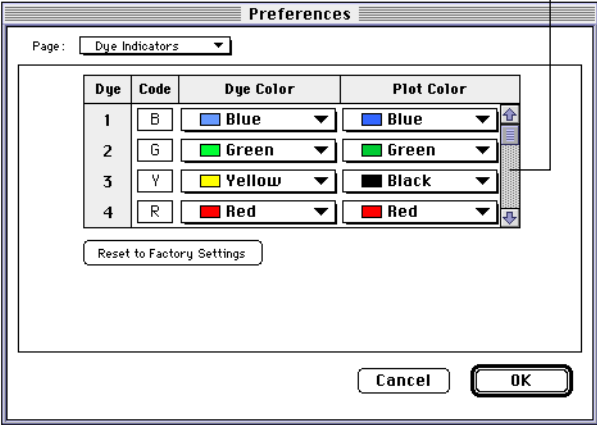
Display or hide by a color by...	Default	Comment
clicking the colored boxes near the top-left corner of the Gel File window. 	All colors displayed.	Any change you make in the button settings is saved in the gel file and used the next time that the file is opened.

Changing Dye Indicators

The following procedure describes how to change the dye indicator colors.

You may want to change the dye colors to make the colors easier to distinguish. For information on changing the default dye indicator colors, see “Setting Dye Indicator Preferences” on page 5-15.

To change the dye indicator colors:

Step	Action					
1	<p>There are two ways to display the Dye Indicators dialog box:</p> <table border="1"> <thead> <tr> <th>You can....</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>Choose Preferences from the Settings menu and Dye Indicators from the submenu.</td> <td rowspan="2">The Dye Indicators dialog box appears.</td> </tr> <tr> <td>Hold the Option key and click one of the colored boxes in the upper corner of the gel display.</td> </tr> </tbody> </table> <p>Note Use the scroll bar to view additional dyes.</p> 	You can....	Result	Choose Preferences from the Settings menu and Dye Indicators from the submenu.	The Dye Indicators dialog box appears.	Hold the Option key and click one of the colored boxes in the upper corner of the gel display.
You can....	Result					
Choose Preferences from the Settings menu and Dye Indicators from the submenu.	The Dye Indicators dialog box appears.					
Hold the Option key and click one of the colored boxes in the upper corner of the gel display.						

To change the dye indicator colors: *(continued)*

Step	Action						
2	Select a color from the pop-up menu to change the Dye color or the Plot color. Changing dye and plot colors has the following affects:						
	<table border="1"> <thead> <tr> <th>Choosing a color from...</th> <th>Shows the colors...</th> </tr> </thead> <tbody> <tr> <td>Dye Color column</td> <td>that represent the dyes in the Analysis Control and the Results Control windows and in the gel image.</td> </tr> <tr> <td>Plot Color column</td> <td>used for displaying the data in electropherograms.</td> </tr> </tbody> </table>	Choosing a color from...	Shows the colors...	Dye Color column	that represent the dyes in the Analysis Control and the Results Control windows and in the gel image.	Plot Color column	used for displaying the data in electropherograms.
	Choosing a color from...	Shows the colors...					
Dye Color column	that represent the dyes in the Analysis Control and the Results Control windows and in the gel image.						
Plot Color column	used for displaying the data in electropherograms.						

Adjusting the Contrast

Use the Adjust Gel Contrast dialog box to increase or reduce the intensity of individual colors in the gel image.

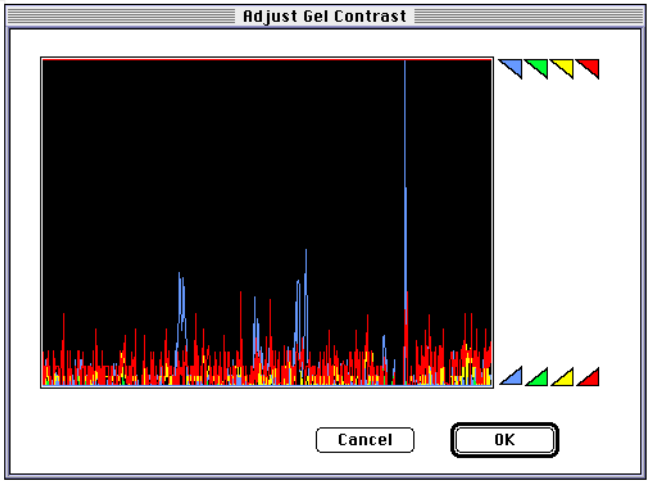
These kinds of adjustments can make it easier to see the data in the gel and can improve the appearance of the gel for presentation.

Note While using the Adjust Gel Contrast dialog box to adjust the contrast, the changes are shown in the background gel.

To adjust the contrast in the gel image:

Step	Action
1	Select a lane in the gel image that contains the color you want to adjust. Note Changes affect the entire gel, not just the selected lane.
2	Choose Adjust Gel Contrast (⌘ J) from the Gel menu. The Adjust Gel Contrast dialog box appears (see below).

To adjust the contrast in the gel image: *(continued)*

Step	Action
	
3	<p>Place the cursor on the triangle for the color you want to adjust, then hold down the mouse button and drag the triangle up or down to a new position.</p> <p>Note The changes take effect in the gel image as you move the triangles.</p>

To adjust the contrast in the gel image: *(continued)*

Step	Action						
<p data-bbox="400 181 419 204">4</p> <p data-bbox="471 181 825 204">You can take the following action.</p> <p data-bbox="471 230 1163 285">Note It is best to adjust one color, apply the change, and view the effect in the gel image before you adjust another color.</p>	<table border="1"> <thead> <tr> <th data-bbox="471 332 825 363">If you want to...</th> <th data-bbox="825 332 1184 363">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="471 363 825 625"> <p data-bbox="481 370 815 393">increase the intensity of a color</p> </td> <td data-bbox="825 363 1184 625"> <p data-bbox="838 370 1130 425">pull the top triangle for that color down.</p> <p data-bbox="838 448 1159 613">For example, pull the top blue triangle down until it is somewhat above the tallest blue data peaks (not primer peaks) in the displayed Slice View.</p> </td> </tr> <tr> <td data-bbox="471 625 825 915"> <p data-bbox="481 631 753 686">suppress the background noise of a color</p> </td> <td data-bbox="825 625 1184 915"> <p data-bbox="838 631 1167 686">pull the bottom triangle for that color up.</p> <p data-bbox="838 709 1167 907">For example, to correct a red background haze because of signal noise or because the signal baseline is not flat, pull the bottom red triangle up until it is just above the baseline and noise in the display.</p> </td> </tr> </tbody> </table>	If you want to...	Then...	<p data-bbox="481 370 815 393">increase the intensity of a color</p>	<p data-bbox="838 370 1130 425">pull the top triangle for that color down.</p> <p data-bbox="838 448 1159 613">For example, pull the top blue triangle down until it is somewhat above the tallest blue data peaks (not primer peaks) in the displayed Slice View.</p>	<p data-bbox="481 631 753 686">suppress the background noise of a color</p>	<p data-bbox="838 631 1167 686">pull the bottom triangle for that color up.</p> <p data-bbox="838 709 1167 907">For example, to correct a red background haze because of signal noise or because the signal baseline is not flat, pull the bottom red triangle up until it is just above the baseline and noise in the display.</p>
	If you want to...	Then...					
	<p data-bbox="481 370 815 393">increase the intensity of a color</p>	<p data-bbox="838 370 1130 425">pull the top triangle for that color down.</p> <p data-bbox="838 448 1159 613">For example, pull the top blue triangle down until it is somewhat above the tallest blue data peaks (not primer peaks) in the displayed Slice View.</p>					
<p data-bbox="481 631 753 686">suppress the background noise of a color</p>	<p data-bbox="838 631 1167 686">pull the bottom triangle for that color up.</p> <p data-bbox="838 709 1167 907">For example, to correct a red background haze because of signal noise or because the signal baseline is not flat, pull the bottom red triangle up until it is just above the baseline and noise in the display.</p>						
<p data-bbox="400 928 419 951">5</p>	<p data-bbox="471 928 1170 984">Click OK to dismiss the dialog box when you are satisfied with the gel contrast.</p>						

Removing Gel Contrast Changes

To remove the changes you can:

- ◆ Choose Undo Adjust Contrast (⌘ Z) from the Edit menu.

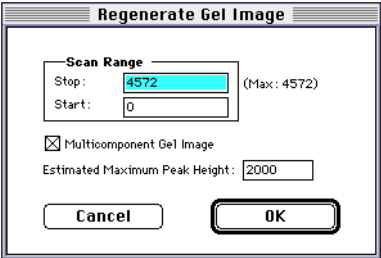
Note Can only undo if you do not choose another command, that is, after adjusting the gel contrast if you choose Undo Adjust Contrast you cannot do segment tracking.

- ◆ Redisplay the Adjust Gel Contrast dialog box and manually reset the settings.
- ◆ Close the Gel File window without saving the changes.

Regenerating the Gel Image

Regenerating the gel image is useful, for example, if there is unusable data near the beginning or the end of the run, the maximum peak height is different from what you expected, or when a new matrix is installed.

To regenerate the gel image:

Step	Action										
1	<p>Choose Regenerate Gel Image from the Gel menu.</p> <p>The Regenerate Gel Image dialog box appears. The values displayed are an example that were used to create the current gel image.</p> 										
2	<p>Make any required changes in the dialog box values, as follows:</p> <table border="1"> <thead> <tr> <th>Use the...</th> <th>To...</th> </tr> </thead> <tbody> <tr> <td>Stop text entry field</td> <td>include the last scan value desired in the gel image. The Max value indicates the total number of scans in the current gel file.</td> </tr> <tr> <td>Start text entry field</td> <td>include the first scan value in the gel image.</td> </tr> <tr> <td>Multicomponent Gel Image checkbox</td> <td>cause GeneScan Analysis Software to apply the assigned matrix file to the gel image. For information on the matrix file, see Chapter 6, "Making a Matrix File."</td> </tr> <tr> <td>Estimated Maximum Peak Height text entry field</td> <td>enter the maximum signal level expected from samples in the run.</td> </tr> </tbody> </table>	Use the...	To...	Stop text entry field	include the last scan value desired in the gel image. The Max value indicates the total number of scans in the current gel file.	Start text entry field	include the first scan value in the gel image.	Multicomponent Gel Image checkbox	cause GeneScan Analysis Software to apply the assigned matrix file to the gel image. For information on the matrix file, see Chapter 6, "Making a Matrix File."	Estimated Maximum Peak Height text entry field	enter the maximum signal level expected from samples in the run.
Use the...	To...										
Stop text entry field	include the last scan value desired in the gel image. The Max value indicates the total number of scans in the current gel file.										
Start text entry field	include the first scan value in the gel image.										
Multicomponent Gel Image checkbox	cause GeneScan Analysis Software to apply the assigned matrix file to the gel image. For information on the matrix file, see Chapter 6, "Making a Matrix File."										
Estimated Maximum Peak Height text entry field	enter the maximum signal level expected from samples in the run.										

To regenerate the gel image: *(continued)*

Step	Action
3	Click OK to close the dialog box and start regenerating the gel image. Note ⌘-(period) cancels the regeneration.

Installing New Matrix Information

About the Matrix File

The Data Collection software copies the matrix information in the specified matrix file to the gel file during data collection. The GeneScan Analysis Software uses this matrix information to multicomponent the gel image (Multicomponenting is the adjusting for spectral overlap of the fluorescent dyes). It also copies this information during lane extraction to each Sample file for use during Sample file analysis.

IMPORTANT If you install new matrix file information after extracting the sample data from the gel file, the Sample files will not contain the new information. You must either regenerate the Sample files, or install the new matrix directly to the existing Sample files.


How to Install a New Matrix File

If you specified the wrong matrix file or no matrix file in data collection, you can install a new matrix file in the gel image. You can also specify a different matrix file in order to test the results of using different matrices on the samples.

To install a new matrix file in a gel image:

Step	Action
1	Open the gel image.
2	Choose Install New Gel Matrix from the Gel menu. A directory dialog box appears. The Folder Preferences settings determine where the GeneScan Analysis Software looks for the matrix file. For more information, see "Specifying File Locations" on page 5-42.

To install a new matrix file in a gel image: *(continued)*

Step	Action
3	<p>Find and select the desired matrix file, then choose Open.</p> <p>The dialog box closes and the information is added to the gel file.</p> <p>If the new matrix installs correctly, the following dialog box appears.</p> <div data-bbox="525 334 1107 594" style="border: 1px solid black; padding: 10px;"><p>Installation of "matix file1" within the Gel File was successful.</p><p>NOTE: Previously extracted sample files DO NOT contain the newly installed matrix.</p><p>Do you want to regenerate the Gel Image?</p><p><input type="button" value="Regenerate Image"/> <input checked="" type="button" value="Don't Regenerate Image"/></p></div>
4	<p>Choose whether or not to regenerate the gel image.</p> <p>Note If you do not regenerate the gel image now, the matrix is included in the data when Sample files are generated.</p> <p>For more information on matrix files, see Chapter 6, "Making a Matrix File."</p>

Verifying Gel Information

Where the Information is Archived

Gel and run information is archived in the following:

- ◆ Data Collection run error log.
- ◆ Gel Info window.
- ◆ Gel image.
- ◆ Gel Sample Sheet.

Data Collection Run Error Log

The Data Collection Run Error log can provide information about Macintosh® computer or instrument errors that occurred during the run.

Gel Info Window

What the Gel Info Window Displays

The window displays information about the run conditions and parameters, and the gel image. You can change some of this information for your record keeping purposes, and other information, such as the date of the run, cannot be changed. The information in this window is saved with each Sample file generated from the gel file.

IMPORTANT Use the Matrix file section in the Gel Info window to verify what matrix is attached to the Gel file and use the Multicomponented section to verify whether or not the Gel file is multicomponented.

How to Display the Gel Info Window

There are two ways to display the Gel Info window:

You can	Result
click the Gel Info button on the Gel File window	The Gel Info window appears (see "Gel Info Window Example" below).
choose Gel Info from the Gel menu	

Gel Info Window Example

Gel file - Gel Info	
Run Information	
User Name: Kathy	Run Date: Tue, Jul 2, 1996
Instrument: 377	Start Time: 7:25:22 AM
Data Coll. Version: 1.1	Run Duration: 2 Hrs 15 Secs
	Total Scans: 4572
Gel Characteristics	
Gel Type:	Number of Channels: 194
Gel Percent: 0.00 %	Gel Thickness: 0.00 mm
Well-To-Read Distance: 36.0 cm	Number of Lanes: 34
	Number of Dyes: 4
Gel Image Information	
Gel Image Range: 0 - 4572	Multicomponented: Yes
Estimated Maximum Peak Height: 2000	Matrix File: matrix file1

Gel Image Review the gel image for a general measure of the quality of the run. For more information, see “How to Adjust the Gel Image” on page 2-18.

IMPORTANT If you make any of the changes to the scan range after extracting the sample data from the gel file, re-extract the Sample files to include the new information. For information, see “Regenerating the Gel Image” on page 2-23.

What to Review in the Gel Image

Review these items...	Take this action...
<ul style="list-style-type: none"> ◆ Inspect the general condition of the bands in the lanes in the gel image. ◆ Are the fluorescent signals displayed as discreet horizontal bands? 	This may indicate a poor gel.
<ul style="list-style-type: none"> ◆ Are any of the colors too bright or too dark? ◆ Is there a green or red haze? ◆ Is this something that you can fix by adjusting the gel image contrast? 	Either: <ul style="list-style-type: none"> ◆ Adjust gel contrast (see page 2-18). ◆ Adjust Estimated Maximum Peak Height (see page 2-6), or ◆ Rerun the gel.
<ul style="list-style-type: none"> ◆ Inspect the lane markers. ◆ Verify that each lane marker corresponds to a sample as designated in the Sample Sheet. 	Adjust the locations of the lane markers, as necessary.

Review these items...	Take this action...
<ul style="list-style-type: none"> ◆ Inspect the tracker lines. ◆ Each tracker line should be centered over the brightest part of the bands it is tracking through. 	Adjust the tracker line placement, as necessary.
<ul style="list-style-type: none"> ◆ Inspect the Slice View for each lane. ◆ Verify that all fragments in the size standard are included in the current tracking. 	Adjust the tracker line placement, as necessary.

Sample Sheet Review the Sample Sheet to confirm that the information in the fields is correct. If necessary, edit the Sample Sheet information (instrument file, sample names, comments, and so forth) that is automatically transferred to the Sample files or to the BioLIMS database.




For more information, see “Displaying the Sample Sheet” on page 2-29.

Displaying the Sample Sheet






What is the Sample Sheet The Sample Sheet contains sample information and is copied to the Data Collection run file before you start the run.

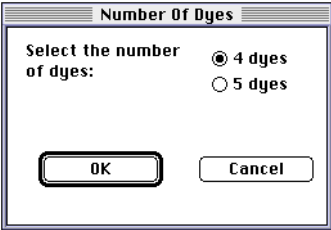
Using the Copy of the Sample Sheet The Sample Sheet copy is automatically copied to the gel file. You can view, edit, and print this Sample Sheet copy whenever the gel file is open. Changes you make in this copy are stored in the gel file; they do not affect the original Sample Sheet file.

Displaying the Copied Sample Sheet To display the copied Sample Sheet:

Step	Action						
1	Open the gel file.						
2	There are two ways to display the Sample Sheet. <table border="1" data-bbox="510 682 1236 860"> <thead> <tr> <th>You can...</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>click the Sample Sheet button </td> <td>The Sample Sheet window appears.</td> </tr> <tr> <td>choose Gel Sample Sheet from the Gel menu</td> <td>See "Sample Sheet Example" on page 2-31.</td> </tr> </tbody> </table>	You can...	Result	click the Sample Sheet button 	The Sample Sheet window appears.	choose Gel Sample Sheet from the Gel menu	See "Sample Sheet Example" on page 2-31.
You can...	Result						
click the Sample Sheet button 	The Sample Sheet window appears.						
choose Gel Sample Sheet from the Gel menu	See "Sample Sheet Example" on page 2-31.						

Displaying a New Sample Sheet To display a new Sample Sheet:

Step	Action
1	Open the GeneScan Analysis Software program.
2	Choose New from the File menu. The Create New dialog box appears. <div data-bbox="524 1193 1095 1356" style="border: 1px solid black; padding: 10px; margin: 10px 0;"> <p>Create New:</p> <div style="display: flex; justify-content: space-around; align-items: center;"> <div style="text-align: center;">  Project </div> <div style="text-align: center;">  Sample Sheet </div> <div style="text-align: center;">  Analysis Parameters </div> <div style="text-align: center;">  Size Standard </div> <div style="text-align: center;">  Matrix </div> <div style="border: 1px solid black; border-radius: 5px; padding: 5px 15px;">Cancel</div> </div> </div>

Step	Action
3	<p>Click the Sample Sheet icon.</p> <p>The following dialog box appears.</p> 
4	<p>Click the radio button that corresponds to the number of dyes contained in your gel file and click OK.</p> <p>An untitled Sample Sheet appears. See "Sample Sheet Example" on page 2-31.</p>

About the Sample Sheet

Sample Sheet Example The following is an example of a Sample Sheet for DNA fragments labeled with a blue dye and the GeneScan-350 Rox Internal Lane Size Standard:

1	2	3	4	5	6	7	8	9	10	11	12
Used	File Name	Sample Name	Dye/Std	Sample Info	Comment	A	P	Collection Name	Collection Comment	Collection Owner	
<input checked="" type="checkbox"/>	sample 1	1347-02	B	1347-02		<input type="checkbox"/>	<input type="checkbox"/>				
			G			<input type="checkbox"/>	<input type="checkbox"/>				
			Y			<input type="checkbox"/>	<input type="checkbox"/>				
			R	Rox 350		<input checked="" type="checkbox"/>	<input type="checkbox"/>				
<input checked="" type="checkbox"/>	sample 2	1347-01	B	1347-01		<input type="checkbox"/>	<input type="checkbox"/>				
			G			<input type="checkbox"/>	<input type="checkbox"/>				
			Y			<input type="checkbox"/>	<input type="checkbox"/>				
			R	Rox 350		<input checked="" type="checkbox"/>	<input type="checkbox"/>				
<input checked="" type="checkbox"/>	sample 3	884-15	B	884-14		<input type="checkbox"/>	<input type="checkbox"/>				
			G			<input type="checkbox"/>	<input type="checkbox"/>				
			Y			<input type="checkbox"/>	<input type="checkbox"/>				
			R	Rox 350		<input checked="" type="checkbox"/>	<input type="checkbox"/>				
<input checked="" type="checkbox"/>	sample 4	884-16	B	884-16		<input type="checkbox"/>	<input type="checkbox"/>				
			G			<input type="checkbox"/>	<input type="checkbox"/>				
			Y			<input type="checkbox"/>	<input type="checkbox"/>				
			R	Rox 350		<input checked="" type="checkbox"/>	<input type="checkbox"/>				
<input checked="" type="checkbox"/>	sample 5	1340-01	B	1340-01		<input type="checkbox"/>	<input type="checkbox"/>				
			G			<input type="checkbox"/>	<input type="checkbox"/>				
			Y			<input type="checkbox"/>	<input type="checkbox"/>				
			R	Rox 350		<input checked="" type="checkbox"/>	<input type="checkbox"/>				
<input checked="" type="checkbox"/>	sample 6	884-17	B	884-17		<input type="checkbox"/>	<input type="checkbox"/>				
			G			<input type="checkbox"/>	<input type="checkbox"/>				
			Y			<input type="checkbox"/>	<input type="checkbox"/>				
			R	Rox 350		<input checked="" type="checkbox"/>	<input type="checkbox"/>				

Sample Sheet Described The following table describes the above diagram.

Sample Sheet described:

No.	Description
1	<p>Lane number for the sample.</p> <p>GeneScan Analysis Software assigns lane numbers to the gel lanes based on the numbers in this column.</p> <p>If a lane in the gel is empty, a corresponding empty row must appear in the Sample Sheet.</p>

Sample Sheet described: *(continued)*

No.	Description						
2	<p>Select this checkbox to mark the corresponding lane in the gel image as Used.</p> <p>When the GeneScan Analysis Software extracts the sample information from the gel file, it creates Sample files only from lanes marked Used.</p> <p>Note Use the Fill Down command from the Edit menu to check all lanes as being "Used".</p> <p>If you change the setting for this checkbox, the corresponding lane setting in the gel file is automatically changed.</p>						
3	<p>Name of the Sample file created for the data in this lane.</p> <p>The Sample file is saved to the hard disk, into the Run folder next to the gel file.</p> <p>If the data is extracted into a BioLIMS database, then this is the name assigned to the database record for that sample file.</p>						
4	<p>Name of the sample, which is used by the Data Collection software to generate the name for the Sample file.</p> <p>The following table lists how the names correspond between the Data Collection software and the GeneScan Analysis Software:</p> <table border="1" data-bbox="467 867 1184 1019"> <thead> <tr> <th data-bbox="467 867 830 906">Data Collection software</th> <th data-bbox="830 867 1184 906">GeneScan Analysis Software</th> </tr> </thead> <tbody> <tr> <td data-bbox="467 906 830 945">Sample Name</td> <td data-bbox="830 906 1184 945">Sample Name</td> </tr> <tr> <td data-bbox="467 945 830 1019">Sample Name and lane number</td> <td data-bbox="830 945 1184 1019">Sample</td> </tr> </tbody> </table> <p>While this information is often the same as the sample info field, it is used by the software for different purposes.</p>	Data Collection software	GeneScan Analysis Software	Sample Name	Sample Name	Sample Name and lane number	Sample
Data Collection software	GeneScan Analysis Software						
Sample Name	Sample Name						
Sample Name and lane number	Sample						
5	<p>Indicates which dye contains the standard.</p>						
6	<p>Information about the sample, usually patient or sample ID.</p> <p>Note This is a required field if the data is to be analyzed using the Genotyper software and stored in GenBase.</p>						
7	<p>Comments about the sample.</p>						
8	<p>Select this checkbox for the GeneScan Analysis Software to automatically analyze the Sample files after generating the gel image.</p>						

Sample Sheet described: *(continued)*

No.	Description
9	When this checkbox and the checkbox labeled A are selected, the GeneScan Analysis Software automatically prints the Sample file information after it automatically analyzes the files.
10	<p>Collection Name—If the samples are extracted into the BioLIMS 2.0 database, this is the name of the collection that will contain them.</p> <p>If a collection of that name does not exist in the database, one is created.</p> <p>See “Editing Collection Information” on page 2-34.</p> <p>Use the Collection Manager to group samples in any number of desired collections (see Appendix E, “Using the BioLIMS Database.”</p>
11	Collection Comment—Comment text associated with BioLIMS 2.0 collection name specified in the Collection Name field described above.
12	Collection Owner—Collection Creator text associated with BioLIMS 2.0 collection name specified in the Collection Name field.

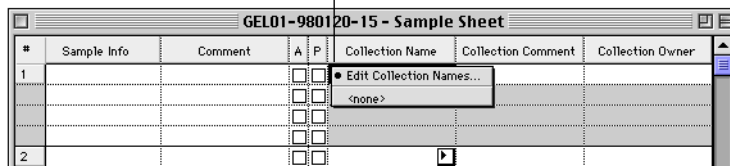
Working with the Sample Sheet

Editing Collection Information

How to Edit Collection Information

Use the pop-up menu in the Collection Name column to edit collection information that corresponds to that sample.

Pop-up menu



Using the Pop-Up Menu

To use the pop-up menu to edit the collection information:

If you want to...	Then...
select a collection	choose a collection from the pop-up menu The Collection Name, Collection comment, and Collection owner fields use this information.
edit a collection	a. Choose the Edit Collection Names pop-up menu. The Collection Name Editor dialog box appears. <div data-bbox="569 945 1166 1344" style="border: 1px solid black; padding: 5px; margin: 10px 0;"> </div> b. Edit the information in the Collection Name, Collection Comment, or Collection Owner fields and click OK.

To use the pop-up menu to edit the collection information: *(continued)*

If you want to...	Then...
add a collection	<ol style="list-style-type: none">a. Choose the Edit Collection Names pop-up menu. The Collection Name Editor dialog box appears (see “edit a collection” above).b. Click Add Row.c. Enter a Collection Name, Collection Comment, and Collection Owner and click OK.
delete a collection	<ol style="list-style-type: none">a. Choose the Edit Collection Names pop-up menu. The Collection Name Editor dialog box appears (see “edit a collection” above).b. Select the row that contains the collection you want to delete.c. Click Delete row and click OK.

Changing Column Widths

To change the width of columns in the Sample Sheet:

Step	Action
1	Place the cursor on the divider line to the right of the column title.
2	When the cursor changes to two arrows, hold down the mouse button and drag the line to the desired location.

Editing the Sample Sheet

To edit the Sample Sheet, you can:

- ◆ Type in new text.

Note Use the Fill Down command to apply a change to all the rows in a column by entering the text in one of the rows, clicking the column header, and then choosing Fill Down from the Edit menu.

- ◆ Use the Edit menu commands.
 - ◆ Select and deselect checkboxes.
-

Printing the Sample Sheet

There are two ways to print the Sample Sheet:

If you choose from the File menu...	Then...
Print (⌘ P)	the standard Print dialog box appears, choose any options, and click Print.
Print One when the Sample Sheet is active	one copy of the selected information is printed, bypassing the standard print dialog box.

Installing a New Sample Sheet

Use the following procedure to replace the original Sample Sheet information by installing a new Sample Sheet.

IMPORTANT If you change the Sample Sheet after extracting Sample files from the gel file, you must re-extract the Sample files to include the new information. Make sure that the checkbox labeled Used is selected for each sample that you want to re-extract.

To install a new Sample Sheet:

Step	Action
1	Choose Install New Sample Sheet from the Gel menu. A directory dialog box appears.
2	Select a new Sample Sheet from the directory dialog box. The new Sample Sheet replaces the original Sample Sheet copied to the gel file.

Importing Data to Sample Sheets

To import data to a Sample Sheet:

Step	Action
1	<p>With the Sample Sheet open, choose Import from the File menu. A directory dialog box appears.</p> <p>Note Ensure that the text file you are importing is complete and properly formatted for a tab-delimited text file, for example a spread sheet program.</p> <p>For more information, refer to the <i>ABI PRISM 377 Instrument User's Manual</i>.</p>
2	<p>Select a text file from the directory dialog box and click Open. The data in the Sample Sheet is replaced by the imported text file.</p>

Exporting Data From Sample Sheets

To export data from Sample Sheets.

Step	Action
1	<p>With the Sample Sheet open, choose Export from the File menu. A directory dialog box appears.</p>
2	<p>Take the following action:</p> <ol style="list-style-type: none">Enter a name for the text file.Choose a folder location to save the file and click Save.

About Displaying Regions of Off-Scale Data

Why Displaying Off-Scale Data is Important

Off-scale data refers to raw fluorescent signals that exceed the linear dynamic range of the CCD camera detection system. A data point is off-scale when the raw fluorescent signal is 8,191 rfu. Peak quantitation of off-scale data is not accurate; this includes both peak area and peak height. Additionally, the matrix file may not be accurately applied, often resulting in pull-up peaks and poor baselining.

Off-Scale Detection and GeneScan

The off-scale detection features of the GeneScan Analysis Software identify and flag such data points in the bands of a gel image and the peaks of an electropherogram. However, these features must be used in conjunction with “suppress left-right averaging” in Data Collection (refer to Setting Preferences, General Settings of chapter 4 in the ABI PRISM 377 User’s Manual) in order to detect all instances of off-scale data.

Displaying Off-Scale Data

Displaying Off-Scale Data In the Gel Image

Take this action	Description
Click the OS button in the Gel File window to display where the data went off-scale data in the gel image. Note To see off-scale portions in the gel image, click the Vertical Expand button. The resolution of the full display is low and does not show the off-scale data.	Each pixel that is off-scale is displayed as a white dot. The other colors in the gel are turned off. When finished viewing the off-scale data, turn on the other colors by clicking each color button or by clicking the OS button again to deselect it.

Displaying Off-Scale Data In the Electropherogram

Pre-Averaging Offscale Detection must be selected when extracting lanes for accurate displays of off-scale data in electropherograms. During the extraction process, each data point used to generate an average value for a scan number (i.e., channel averaging) is compared to the off-scale value of 8191.

Note If you extracted a sample file with the Pre-Averaging Offscale Detection checkbox selected, then pre-averaging has a value of zero in the Sample Info View.

How Pre-Averaging Offscale Detection Works

The following table lists how the Pre-Averaging Offscale Detection feature identifies off-scale data points:

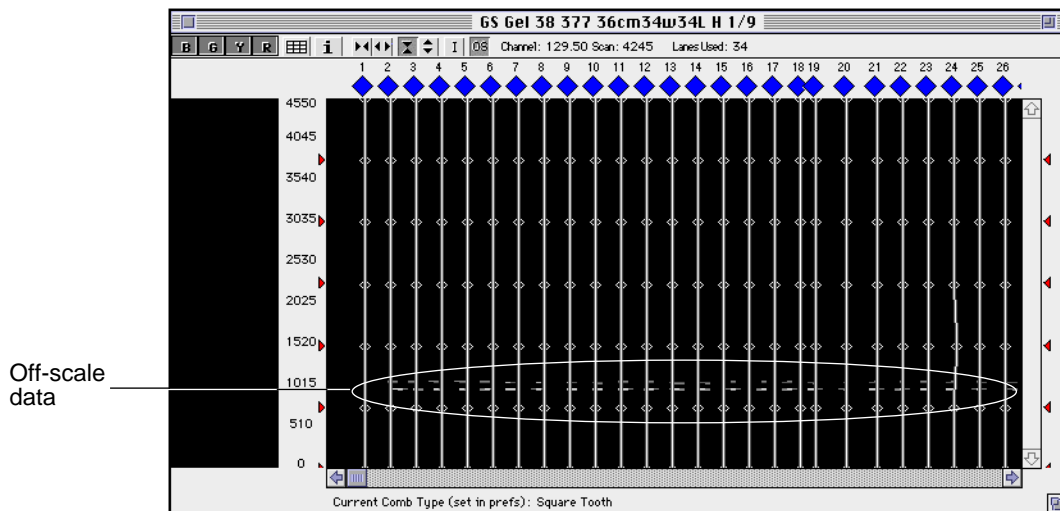
If...	Then...
each data point used in averaging is less than 8191 rfu (relative fluorescent units)	the peak in the electropherogram is not off-scale and appears normal.
any one of the data points used in the averaging is 8191 rfu	<p>no averaging takes place and 8191 rfu is written for that scan number of that lane.</p> <p>In the electropherogram, the off-scale regions are displayed by drawing red vertical regions where the off-scale data is present.</p> <p>How to Display the Red Regions</p> <ol style="list-style-type: none">Choose Preferences from the Settings menu and Results Display from the submenu.Select the checkbox labeled "Show Offscale Regions." <p>Note You can choose Hide/Show Offscale Regions from the View menu if the checkbox labeled Show Offscale Regions is not selected.</p>

About Flat Topped Peaks

An additional feature is peaks that contain off-scale data points are drawn in the electropherograms as "flat topped." That is, the top section of the peak is flat rather than pointed (see "Electropherogram Displaying the Flat Topped Effect" on page 2-42).

This feature may be seen when the data is analyzed with no or light smoothing; the flat topped peaks may not be apparent with heavy smoothing. Choose Analysis Parameters from the Settings Menu (see "Smooth Options" on page 5-20).

Gel File Window Example The following is an example of off-scale data displayed in a Gel File window. Each white line contains off-scale data points.

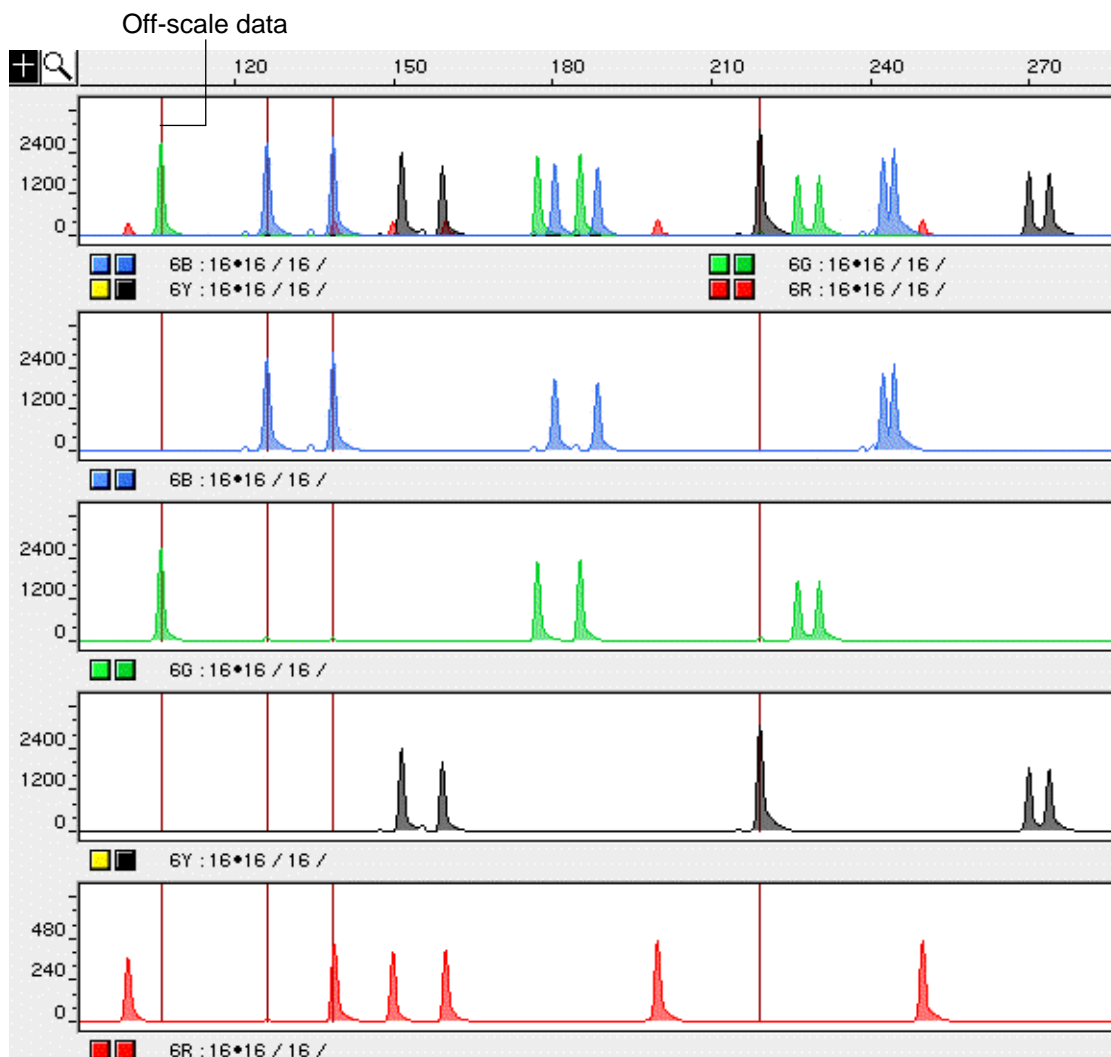


Electropherogram Examples

The following are two examples of off-scale data displayed in electropherograms. Each red bar/region indicates an off-scale data point.

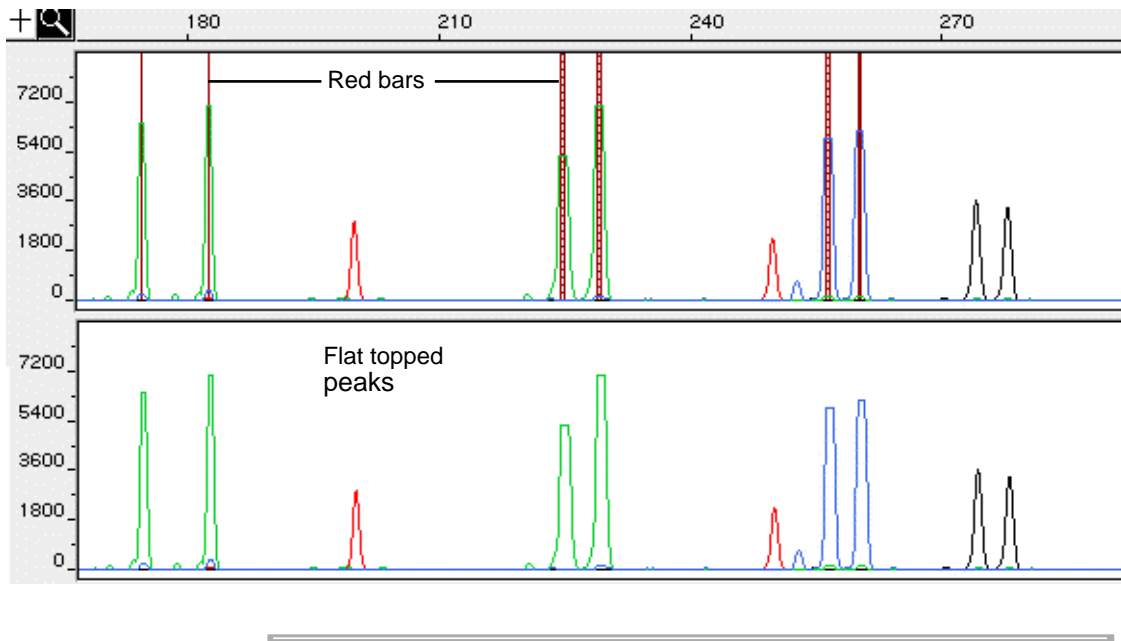
Electropherogram Displaying Off-Scale Data

Note Choose the Zoom In command from the View menu to more clearly show the areas of off-scale data.



Electropherogram Displaying the Flat Topped Effect

The off-scale peaks in the zoomed-in electropherograms illustrate the flat topped effect. The second electropherogram has the red bars removed.



Working with Lane Markers

Introduction When the GeneScan Analysis Software first opens a gel file, it adds lane numbers, lane markers, and tracker lines to the gel image. In most cases, if you used a good gel for your run and the run settings are correct, the gel file should be properly tracked. If, however, the tracking is incorrect because the signal is weak and the tracker missed a lane, you might need to make some changes.

Topics in This Section This section includes the following topics:

For this topic	See page
Moving Between Lanes and Channels	2-44
Using Control Points	2-45
Adding and Deleting Control Rows	2-46
What Marking Lanes as Used/Unused Means	2-47
Marking Lanes as Used or Unused	2-47
Moving Mislaced Lane Markers	2-48
Lane Marker Rules	2-50
About Marking Lanes for Extraction	2-50
Marking All Lanes for Extraction	2-51
Unmarking All Lanes for Extraction	2-51
Marking a Single Lane for Extraction	2-52
Unmarking a Single Lane for Extraction	2-52

Moving Between Lanes and Channels

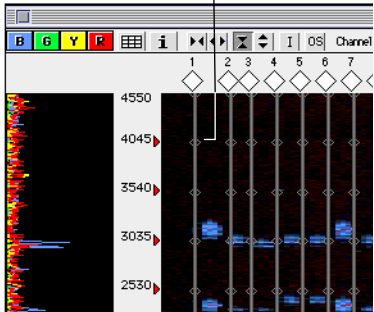
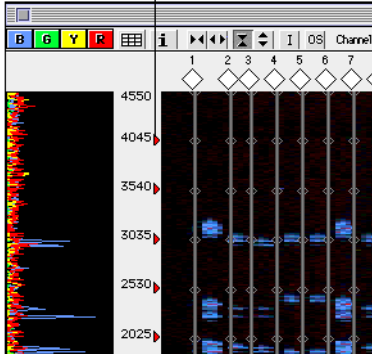
The following table lists how to move between lanes and channels:

If you want to move...	Then...									
quickly from one lane to the next lane	<p>You can take the following action:</p> <table border="1" data-bbox="706 293 1184 571"> <thead> <tr> <th data-bbox="706 293 951 334">To move...</th> <th data-bbox="956 293 1184 334">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="706 337 951 431">the lane</td> <td data-bbox="956 337 1184 431">click the Lane marker of the lane you want to move.</td> </tr> <tr> <td data-bbox="706 435 951 500">one lane to the right</td> <td data-bbox="956 435 1184 500">press the Tab key.</td> </tr> <tr> <td data-bbox="706 503 951 571">one lane to the left</td> <td data-bbox="956 503 1184 571">press the Shift-Tab key.</td> </tr> </tbody> </table>		To move...	Then...	the lane	click the Lane marker of the lane you want to move.	one lane to the right	press the Tab key.	one lane to the left	press the Shift-Tab key.
To move...	Then...									
the lane	click the Lane marker of the lane you want to move.									
one lane to the right	press the Tab key.									
one lane to the left	press the Shift-Tab key.									
the location of a selected lane by one channel	<p>You can take the following action:</p> <ul style="list-style-type: none"> ◆ Click the control point and drag, or ◆ Use the Arrow keys. <p>Note Pressing the Option key while using the Arrow keys moves the control point in 0.1 channel increments.</p>									
a lane segment	Use the arrows to move the segment.									
from one tracker line segment to the next segment	<p>You can take the following action:</p> <table border="1" data-bbox="706 927 1184 1073"> <thead> <tr> <th data-bbox="706 927 951 967">To move...</th> <th data-bbox="956 927 1184 967">Use the...</th> </tr> </thead> <tbody> <tr> <td data-bbox="706 971 951 1003">up one segment</td> <td data-bbox="956 971 1184 1003">Up Arrow key.</td> </tr> <tr> <td data-bbox="706 1006 951 1073">down one line segment</td> <td data-bbox="956 1006 1184 1073">Down Arrow key.</td> </tr> </tbody> </table>		To move...	Use the...	up one segment	Up Arrow key.	down one line segment	Down Arrow key.		
To move...	Use the...									
up one segment	Up Arrow key.									
down one line segment	Down Arrow key.									

Using Control Points

When a gel is tracked, the tracker places a tracker line with spline control points over the strongest signals for its lane. Use the control points to control shape and position the splines. When a control point is selected it appears as a filled in square.

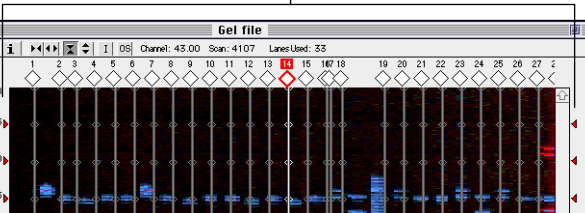
The following table lists how you can use the control points:

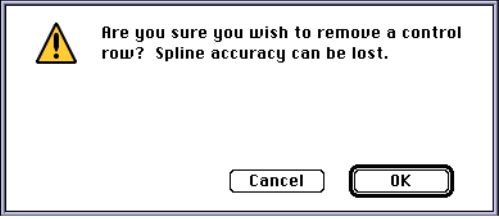
If you want to...	Then...
<p>move a single control point</p> <p style="text-align: center;">Control point</p> 	<ul style="list-style-type: none"> ◆ Click the control point and drag, or ◆ Use the Arrow keys. <p>Note Pressing the Option key while using the Arrow keys moves the control point in 0.1 channel increments.</p>
<p>select an entire row</p> <p style="text-align: center;">Red triangle</p> 	<p>click one of the red triangles at the sides of the gel view.</p> <p>Use the Right or Left Arrow keys to move all the control points in that direction.</p>
<p>select multiple control points</p>	<p>draw a box around a set of control points.</p> <p>All the control points that you drew a box around become enabled.</p> <p>Use the Arrow keys to move all the control points.</p>

If you want to...	Then...
move up or down the control points on the spline	use the Arrow keys.
move forward (right) a lane Note When moving forward the control point is deselected.	press the Tab key.
move left (backwards) a lane Note When moving backwards the control point is deselected.	press the Shift-Tab key.

Adding and Deleting Control Rows

To add and delete control rows:

If you want to...	Then...
add a control row	<p>a. Place the arrow cursor outside the image (left or right side)</p> <p>Place arrow cursor on left or right side</p>  <p>The cursor changes to a symbol and a horizontal line will be displayed across the gel image.</p> <p>b. Click the mouse button to add a control row at the desired location.</p>

If you want to...	Then...
delete a control row	<p>a. Option-click once on the control row triangle that you want to delete.</p> <p>The following dialog box appears.</p> <div data-bbox="669 272 1166 488" style="border: 1px solid black; padding: 10px; margin: 10px auto; width: fit-content;">  <p style="text-align: center;">Are you sure you wish to remove a control row? Spline accuracy can be lost.</p> <p style="text-align: right;"> <input type="button" value="Cancel"/> <input type="button" value="OK"/> </p> </div> <p>b. Click OK.</p>

What Marking Lanes as Used/Unused Means

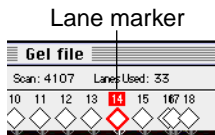
The GeneScan Analysis Software only tracks a lane and extracts sample data if the lane is marked Used in the Gel Sample Sheet. When a lane is marked Used, its lane marker (◆) is colored blue, white, or yellow. All unused lanes have gray lane markers.

By marking lanes as Used or Unused, you specify which lanes contain sample data. This allows the software to correctly number the Used lanes and put the extracted sample data from each lane into the correct Sample file. It also ensures that Sample files are generated from only the Used lanes, and not from empty lanes.

Marking Lanes as Used or Unused

Note Using any of these three methods to mark a lane used or unused changes the setting in both the gel image and the Sample Sheet.

The three ways to mark a lane as used or unused are by:

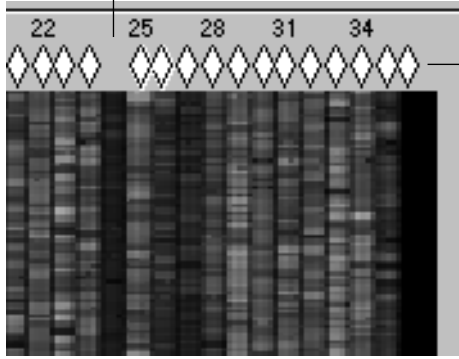


- ◆ Clicking the lane marker, then choosing Mark Lane Used/Unused from the Gel menu.
- ◆ Holding down the ⌘ key while clicking the lane marker.
- ◆ Clicking the Sample Sheet button to open the Sample Sheet that is attached to the gel file, and then selecting or deselecting the Used checkbox for the lane.

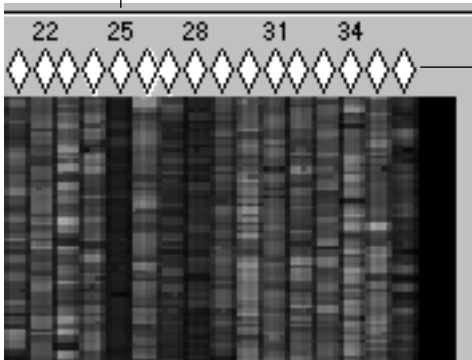
Moving Misplaced Lane Markers

The following procedure describes how to move misplaced lane markers so that the lane numbers are properly aligned with the actual sample lanes. The lane data is written to the correct files with the Sample files are regenerated.

To move misplaced lane markers:

Step	Action
1	<p>Inspect the gel image for incorrectly labeled lanes.</p> <p>For example, in the following illustration, the GeneScan Analysis Software missed the faint signals from lane #25. As a result, lanes #26–36 are mislabeled, and the lane #36 marker is over an unused area to the right of the lanes.</p>  <p>Missed lane #25 has no marker</p> <p>Lane #36 marker is not over a lane</p>

To move misplaced lane markers: *(continued)*

Step	Action
2	<p data-bbox="517 178 1223 267">Click the incorrectly placed marker to select it (the selected marker has a red border), and while holding down the mouse button, drag the lane marker to the correct location.</p> <p data-bbox="517 284 1081 316">All the affected markers are renumbered accordingly.</p> <p data-bbox="517 332 1202 414">Lane markers always remain in numerical order from left to right and are attached to their respective tracker line within a few channels to either side.</p> <p data-bbox="517 430 1216 519">For example, if you drag the #36 marker (in the preceding illustration) to the real lane 25, lanes #25–36 all become correctly marked (as shown in the figure below).</p> <div data-bbox="517 544 1216 966"><p data-bbox="1014 552 1176 600">Lane #25 now has a marker</p><p data-bbox="1014 625 1209 722">Lane #36 marker is now above right-most lane of gel</p></div>

Marking and Unmarking Lanes for Extraction

Lane Marker Rules The GeneScan Analysis Software marks lanes according to the following rules:

If...	Then...	Using...
the lane is identified with sample names in the sample sheet of the Data Collection software	it is automatically marked as Used.	blue markers
the lane is unidentified	it is marked as Unused.	gray markers
you are opening a gel file for the first time	all Used lanes are marked for extraction.	white markers
you have extracted the lanes	each lane is unmarked for extraction.	blue markers
you have edited a lane	it is automatically marked for extraction.	white markers
the lane was inferred by the Tracker software.	the lane is marked.	orange markers
a lane is modified and then extracted and its edited tracking information is not saved to a gel file	this serves as a warning that the latest generated sample does not reflect the saved tracking information.	yellow markers

About Marking Lanes for Extraction During the extraction process, GeneScan Analysis Software only extracts data from gel lanes that are marked for extraction. This allows you to control which lanes to extract when using the Extract Lanes command.

For more information, see “Extracting Data Without Changing the Current Tracker” on page 2-64.

Marking All Lanes for Extraction

To mark all lanes for extraction:

From the Gel menu choose...	Result
Mark All Lanes for Extraction	<p>The color of all markers for Used lanes changes to white.</p> <p>If you choose the Extract Lanes command, the GeneScan Analysis Software uses the current tracker line location to extract the data in all Used lanes and puts the extracted data in new sample files.</p>

Unmarking All Lanes for Extraction

To unmark all lanes for extraction:

Step	Action						
1	Select Unmark All Lanes For Extraction from the Gel menu. The color of all the lane markers changes to blue.						
2	You can take the following action. <table border="1" data-bbox="512 805 1239 989"> <thead> <tr> <th>To...</th> <th>See...</th> </tr> </thead> <tbody> <tr> <td>extract all lanes marked as Used</td> <td>step 2 on page 2-65.</td> </tr> <tr> <td>remark selected lanes for extraction</td> <td>"Marking a Single Lane for Extraction" on page 2-52.</td> </tr> </tbody> </table>	To...	See...	extract all lanes marked as Used	step 2 on page 2-65.	remark selected lanes for extraction	"Marking a Single Lane for Extraction" on page 2-52.
To...	See...						
extract all lanes marked as Used	step 2 on page 2-65.						
remark selected lanes for extraction	"Marking a Single Lane for Extraction" on page 2-52.						

Marking a Single Lane for Extraction

To mark only selected lanes or to extract a lane that was not automatically extracted:

Step	Action						
1	Click the lane marker you want to mark for extraction.						
2	There are two ways to mark a single lane for extraction: <table border="1" data-bbox="467 370 1190 727"> <thead> <tr> <th>You can either...</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>choose Mark Lane for Extraction from the Gel menu</td> <td>The color of the lane marker changes to white.</td> </tr> <tr> <td>press the Option key and click the lane marker</td> <td>When you choose the Extract Lanes command, the GeneScan Analysis Software uses the current tracker line locations to extract the data in this and other similarly marked files, and puts the extracted data into new Sample files.</td> </tr> </tbody> </table>	You can either...	Result	choose Mark Lane for Extraction from the Gel menu	The color of the lane marker changes to white.	press the Option key and click the lane marker	When you choose the Extract Lanes command, the GeneScan Analysis Software uses the current tracker line locations to extract the data in this and other similarly marked files, and puts the extracted data into new Sample files.
You can either...	Result						
choose Mark Lane for Extraction from the Gel menu	The color of the lane marker changes to white.						
press the Option key and click the lane marker	When you choose the Extract Lanes command, the GeneScan Analysis Software uses the current tracker line locations to extract the data in this and other similarly marked files, and puts the extracted data into new Sample files.						

Unmarking a Single Lane for Extraction

To unmark a lane that is marked for extraction:

Step	Action					
1	Click the lane marker for the lane that you want to unmark.					
2	There are two ways to unmark a single lane for extraction: <table border="1" data-bbox="467 1016 1190 1193"> <thead> <tr> <th>You can...</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>choose Unmark Lane for Extraction from the Gel menu</td> <td rowspan="2">The color of the lane marker changes to blue.</td> </tr> <tr> <td>press the Option key and click the lane marker</td> </tr> </tbody> </table>	You can...	Result	choose Unmark Lane for Extraction from the Gel menu	The color of the lane marker changes to blue.	press the Option key and click the lane marker
You can...	Result					
choose Unmark Lane for Extraction from the Gel menu	The color of the lane marker changes to blue.					
press the Option key and click the lane marker						

Working with Tracker Lines

Topics in This Section This section includes the following topics:

Topics	See page
Hiding/Showing Tracker Lines	2-53
Why Positioning Each Lane is Important	2-53
Causes to Misinterpreting Lane Positions	2-54
Verifying Optimal Channel Tracking	2-55
Reviewing Tracker Lines	2-55
Moving Tracker Lines	2-56
Deselecting a Lane	2-56

Hiding/Showing Tracker Lines The following table lists how to hide and show tracker lines. The first time the GeneScan Analysis Software opens a gel file, tracker lines are added to the gel image.

To...	Then...						
turn off the tracker lines	choose Hide Tracker Lines from the Gel Menu. All unselected tracker lines disappear. If a lane is selected, the white tracker line for that lane remains visible.						
turn on the tracker lines	choose Show Tracker Lines from the Gel menu. The following table lists how all the tracker lines are displayed.						
	<table border="1"> <thead> <tr> <th>If...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>no lane is selected</td> <td>all lanes are grey.</td> </tr> <tr> <td>a lane is selected</td> <td>that lane is white and all the other lanes are grey.</td> </tr> </tbody> </table>	If...	Then...	no lane is selected	all lanes are grey.	a lane is selected	that lane is white and all the other lanes are grey.
If...	Then...						
no lane is selected	all lanes are grey.						
a lane is selected	that lane is white and all the other lanes are grey.						
display one tracker line	click the lane marker (◆) for the tracker line you want to view. The unselected tracker lines are greyed out.						

**Why Positioning
Each Lane is
Important**

The GeneScan Analysis Software normally calculates the data values for each lane by averaging the data from selected channels.

Unless each lane is positioned at the optimum channel of the fragment band, the resulting calculations for peak height and peak area may be incorrect. Size calling, however, is still correct.

**Causes to
Misinterpreting
Lane Positions**

The following table lists the possible conditions that may cause automatic tracking to misinterpret lane positions or fail to follow the path of a lane completely.

Condition	Result
Weak signals	<ul style="list-style-type: none">◆ Causes the software to completely miss or be unable to follow a lane, especially if the gel ran abnormally.◆ Although the software creates a track for each Used lane on the gel, the tracker lines might be incorrectly placed, indicating lane positions that do not exist or that are located elsewhere.
Gels that contain extremely bent lanes	<ul style="list-style-type: none">◆ Causes the software to lose track of a lane at some point along the lane's path.◆ Make identification of lane positions difficult.
Failure to complete the sample sheet	<ul style="list-style-type: none">◆ The software uses the Used lane information to determine if what is found corresponds to what was loaded according to the Sample Sheet.◆ If too few or too many lanes are marked Used, then the software must guess at which lanes to throw out and which to keep.
Wrong or bad matrix	The software needs a proper matrix to determine the centers of the bands of the gel.

Verifying Optimal Channel Tracking

To verify optimal channel tracking, examine peak heights in the Slice View of the Gel File window. These peaks represent the selected lane and show the peak intensity.

If the lane is...	Then...
correctly tracked	the white tracker line should appear in the middle of each band.
optimally located	you may need to adjust the line and re-extract the affected lane. For more information, see "Tracking and Extracting Data" on page 2-61.

Reviewing Tracker Lines

To review tracker line placement:

Step	Action
1	In the Gel File window, click the lane marker for the lane you want to review. Note From the currently selected lane, you can also press the Tab key to move right, or press Shift-Tab to move left.
2	Inspect the placement of the tracker line on the lane. <ul style="list-style-type: none">◆ Does the tracker line seem to follow the brightest part of the lane (usually the center) for the full length of the lane?◆ Does the lane drift to the side while the tracker line does not?
3	Inspect the Slice View of the lane data. <ul style="list-style-type: none">◆ Are the raw data peaks high enough for the full length of the lane?◆ Are some peaks very low because the tracker line is not located correctly on the lane?

Moving Tracker Lines

Tracker lines can be moved in the following ways:

To move...	Then...
the line over one channel	a. Click the lane marker for the tracker line you want to move. The lane marker becomes outlined in red, and the tracker lines for all other lanes become hidden or grayed out. b. Press the Left Arrow or Right Arrow key to move the line over one channel.
selected lines to the right	choose Force Selected Lanes to the Right from the Gel menu.
the line in channel increments of 0.1	a. Click the lane marker for the tracker line you want to move. b. Press the Option key and either the Left or Right Arrow.

Deselecting a Lane


To de-select an edited lane:

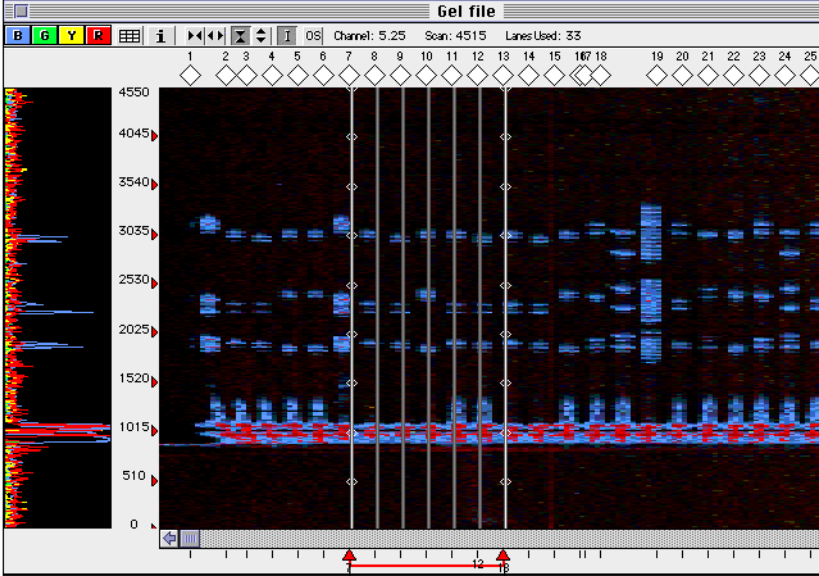
In the lane marker area at the top of the window....	Result						
click the background (between the diamonds that mark the lanes) to deselect the lane marker.	<table border="1"> <thead> <tr> <th>If...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>all tracker lines are displayed</td> <td>all lines become grey.</td> </tr> <tr> <td>all lines are hidden</td> <td>the deselected line becomes hidden.</td> </tr> </tbody> </table>	If...	Then...	all tracker lines are displayed	all lines become grey.	all lines are hidden	the deselected line becomes hidden.
	If...	Then...					
	all tracker lines are displayed	all lines become grey.					
all lines are hidden	the deselected line becomes hidden.						

Interpolating Tracker Lines

Why Use Interpolation Mode Use the interpolation mode to adjust the tracking lines between two vertical points in the gel image. This mode is very useful to manually manipulate the tracker lines for a set of lanes than have anomalous but similar migration patterns. Once the two points are chosen, you can use the control points on the tracker lines to move the lines.

Procedure To use the interpolation mode:

Step	Action
1	With the Gel File window displayed, press the Interpolation button to enter interpolation mode. 
2	In the region being tracked, click the left-hand lane marker at the bottom of the tracker line and then click the right-hand lane marker. These two lanes are the interpolation guides and the lanes between them are interpolated. Note While in interpolation mode, all the other lanes are hidden and no editing of the tracker lines is allowed.
3	Press the Interpolation button again to exit the interpolation mode.

Step	Action						
	 <p style="text-align: center;">Interpolation guides</p>						
4	<p>You can take the following action:</p> <table border="1" data-bbox="467 824 1185 1386"> <thead> <tr> <th data-bbox="467 824 830 863">To...</th> <th data-bbox="830 824 1185 863">You can...</th> </tr> </thead> <tbody> <tr> <td data-bbox="467 863 830 1221">use the control points to move the tracker lanes</td> <td data-bbox="830 863 1185 1221"> <ul style="list-style-type: none"> ◆ Click the control point and drag, or ◆ Use the Arrow keys. <p>Note Pressing the Option key while using the Arrow keys moves the control point in 0.1 channel increments.</p> <p>For more information, see “Using Control Points” on page 2-45.</p> </td> </tr> <tr> <td data-bbox="467 1221 830 1386">track and extract lanes</td> <td data-bbox="830 1221 1185 1386"> <p>choose Extract Lanes from the Gel menu.</p> <p>For more information, see “Tracking Lanes and Extracting Data” on page 2-59.</p> </td> </tr> </tbody> </table>	To...	You can...	use the control points to move the tracker lanes	<ul style="list-style-type: none"> ◆ Click the control point and drag, or ◆ Use the Arrow keys. <p>Note Pressing the Option key while using the Arrow keys moves the control point in 0.1 channel increments.</p> <p>For more information, see “Using Control Points” on page 2-45.</p>	track and extract lanes	<p>choose Extract Lanes from the Gel menu.</p> <p>For more information, see “Tracking Lanes and Extracting Data” on page 2-59.</p>
To...	You can...						
use the control points to move the tracker lanes	<ul style="list-style-type: none"> ◆ Click the control point and drag, or ◆ Use the Arrow keys. <p>Note Pressing the Option key while using the Arrow keys moves the control point in 0.1 channel increments.</p> <p>For more information, see “Using Control Points” on page 2-45.</p>						
track and extract lanes	<p>choose Extract Lanes from the Gel menu.</p> <p>For more information, see “Tracking Lanes and Extracting Data” on page 2-59.</p>						

Tracking Lanes and Extracting Data

Ways to Track and Extract Data **Note** The only action that changes the tracking is choosing the correct comb type (see “How to Set Gel Processing Preferences” on page 2-3). Changes to the Sample Sheet do not effect the tracking.

There are three ways to track and extract data in the gel file:

To...	Choose from the Gel Menu	Comment	See...
track the gel without extracting Sample file data	Track Lanes command to track the gel file based on the current Sample Sheet information.	Both these procedures generate the same tracking lines.	“Tracking the Gel Without Extracting Data” on page 2-60.
both track the lanes and extract data into the Sample files or the BioLIMS database records	Track and Extract Lanes command to calculate the tracking and then extract the Sample file from the tracked gel.		“Tracking and Extracting Data” on page 2-61.
extract Sample file data without changing the current line tracking	Extract Lane command each time you change the tracker line positioning in a lane, or the information in the Sample Sheet, to incorporate the new information.		“Extracting Data Without Changing the Current Tracker” on page 2-64.

If the Run Folder is Not Specified Each time the GeneScan Analysis Software tracks and extracts a gel, the software looks for the data collection Run folder that is specified in the gel file.

If the...	Then...
<ul style="list-style-type: none"> ◆ Software does not locate the folder, or if the 	the program creates a new one in the folder that contains the gel file.
<ul style="list-style-type: none"> ◆ Gel file does not reference one 	The program places new Sample files in the Run folder.

Tracking the Gel Without Extracting Data

Why Use this Procedure

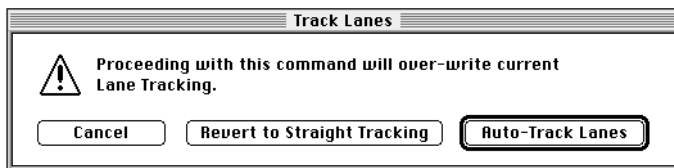
Use this procedure to:

- ◆ View the results of auto-tracking and, if necessary, to correct any tracking errors or tracker line interpolation, before the software extracts the lanes and generates Sample files.
- ◆ Edit tracker information, and to redraw the lines based on original tracking information.

Procedure

Choose Track Lanes from the Gel menu. The Track Lanes dialog box appears.

Note Pressing ⌘-period cancels tracking.



You can take the following actions:

If you choose...	Then this...
Cancel	cancels the tracking operation if you do not want to lose the current tracking information.
Revert to Straight Tracking	adds straight, evenly spaced tracker lines to the gel. For information about moving these lines to the region of the highest signal within the lane, see "Working with Tracker Lines" on page 2-53.

If you choose...	Then this...
<p>Auto-Track Lane</p> <p>IMPORTANT Before choosing Auto-Track Lane, verify that the comb type is set correctly by choosing Preferences from the Settings menu and Gel Preferences from the submenu.</p>	<p>places a tracker line with spline control points over the strongest signals for its lane.</p> <p>For information about:</p> <ul style="list-style-type: none"> ◆ Setting the lane assignment confidence value, see “Confidence Threshold text entry field” on page 2-10. ◆ Moving these lines to the region of the highest signal within the lane, see “Working with Tracker Lines” on page 2-53.

Tracking and Extracting Data

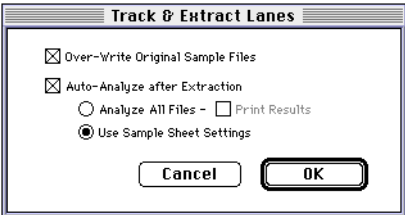
Why Use This Procedure

Use this procedure to automatically track and extract lanes.

IMPORTANT When you Track and Extract a gel file, ensure the Sample Sheet associated with the gel file has the checkbox labeled Used selected for each sample you want to extract. The software only extracts data from used lanes (see “Sample Sheet Example” on page 2-31).

Procedure


To track and extract the gel file:

Step	Action
1	<p>Choose Track & Extract Lanes from the Gel menu.</p> <p>The Track & Extract Lanes dialog box appears.</p> 

To track and extract the gel file: *(continued)*

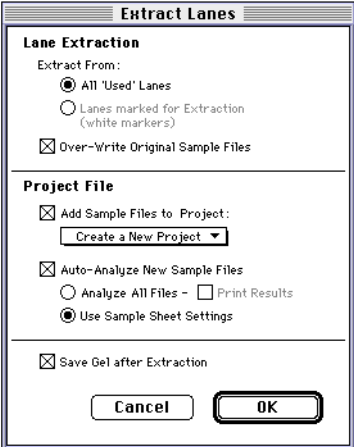
Step	Action						
2	<p data-bbox="471 183 1170 207">Use the Over-Write Original Sample Files checkbox as follows:</p> <table border="1" data-bbox="467 250 1188 646"> <thead> <tr> <th data-bbox="471 256 612 289">If you...</th> <th data-bbox="612 256 1184 289">Then this...</th> </tr> </thead> <tbody> <tr> <td data-bbox="471 297 612 548">select the checkbox</td> <td data-bbox="612 297 1184 548"> <p data-bbox="626 297 1170 354">overwrites any existing Sample files that have the same name with the same sample data.</p> <p data-bbox="626 370 1170 540">Tracking and extracting the gel file a second time without overwriting the original Sample files, creates new Sample files with a dot and a number (.#) appended to the original names. The number increments each time the program creates new files.</p> </td> </tr> <tr> <td data-bbox="471 557 612 638">do not select the checkbox</td> <td data-bbox="612 557 1184 638"> <p data-bbox="626 557 1170 613">preserves Sample file names that already exist and create new Sample files for this data.</p> </td> </tr> </tbody> </table> <p data-bbox="471 686 1170 768">Note In BioLIMS mode, the file name must be unique. If the file name is not unique, then the sample file will not be written to the database.</p> <p data-bbox="471 792 1170 963">Note If a series of numbered files exist, and you discard one of the files and then have the GeneScan Analysis Software re-extract that sample, the software adds the first available number to the new file. For example, if SAMPLE.1, SAMPLE.2, and SAMPLE.3 exist, and you discard SAMPLE.2, GeneScan names the next file SAMPLE.2 (not SAMPLE.4).</p>	If you...	Then this...	select the checkbox	<p data-bbox="626 297 1170 354">overwrites any existing Sample files that have the same name with the same sample data.</p> <p data-bbox="626 370 1170 540">Tracking and extracting the gel file a second time without overwriting the original Sample files, creates new Sample files with a dot and a number (.#) appended to the original names. The number increments each time the program creates new files.</p>	do not select the checkbox	<p data-bbox="626 557 1170 613">preserves Sample file names that already exist and create new Sample files for this data.</p>
If you...	Then this...						
select the checkbox	<p data-bbox="626 297 1170 354">overwrites any existing Sample files that have the same name with the same sample data.</p> <p data-bbox="626 370 1170 540">Tracking and extracting the gel file a second time without overwriting the original Sample files, creates new Sample files with a dot and a number (.#) appended to the original names. The number increments each time the program creates new files.</p>						
do not select the checkbox	<p data-bbox="626 557 1170 613">preserves Sample file names that already exist and create new Sample files for this data.</p>						
3	<p data-bbox="471 979 1170 1036">Use the Auto-Analyze after Extract checkbox to Analyze the new Sample files after extraction is finished.</p> <p data-bbox="471 1052 1170 1109">Do not select this checkbox if you want the Sample files extracted but not analyzed.</p>						
4	<p data-bbox="471 1125 1170 1174">Use the Analyze All Files radio button to analyze all the new Sample files created from the gel file.</p>						
5	<p data-bbox="471 1190 1170 1247">Use the Print Results checkbox to print the results for all new Sample files after analysis.</p>						
6	<p data-bbox="471 1263 1170 1344">Choose Use Sample Sheet Settings radio button to analyze and print only those files that are marked for analysis and printing in the Sample Sheet for the gel file.</p> <p data-bbox="471 1369 1170 1417">Note Select this option if you want to process only some of the new Sample files created from the gel file.</p>						

To track and extract the gel file: *(continued)*

Step	Action								
7	Click OK. See the table below for the action you can take.								
If a project is...	Then...								
open	<p>The following dialog box appears.</p>  <p>And you can take the following action:</p> <table border="1"> <thead> <tr> <th>To...</th> <th>Click...</th> </tr> </thead> <tbody> <tr> <td>add the new Sample files to the open project</td> <td>Use Open Project.</td> </tr> <tr> <td>close the project and create one for the new Sample files</td> <td>Close it.</td> </tr> <tr> <td>cancel the track and extract command</td> <td>Cancel.</td> </tr> </tbody> </table>	To...	Click...	add the new Sample files to the open project	Use Open Project.	close the project and create one for the new Sample files	Close it.	cancel the track and extract command	Cancel.
To...	Click...								
add the new Sample files to the open project	Use Open Project.								
close the project and create one for the new Sample files	Close it.								
cancel the track and extract command	Cancel.								
not open	The GeneScan Analysis Software creates a new project and adds the new Sample files to it.								

Extracting Data Without Changing the Current Tracker

To extract the data without changing current tracker lines:

Step	Action
1	<p>Choose Extract Lanes (⌘ L) from the Gel menu and the Extract Lanes dialog box appears.</p> <p>Note Press ⌘-period (.) to cancel the Extract Lanes operation.</p> 

To extract the data without changing current tracker lines: *(continued)*

Step	Action								
2	In the Lane Extraction portion of the dialog box you can take the following action:								
	<table border="1"> <thead> <tr> <th data-bbox="510 289 880 321">Choose...</th> <th data-bbox="880 289 1239 321">To...</th> </tr> </thead> <tbody> <tr> <td data-bbox="510 321 880 391">All Used Lanes radio button</td> <td data-bbox="880 321 1239 391">generate a new Sample file for every lane marked Used.</td> </tr> <tr> <td data-bbox="510 391 880 488">Lanes marked for Extraction radio button (white markers)</td> <td data-bbox="880 391 1239 488">generate a new Sample file only for each lane with a white lane marker.</td> </tr> <tr> <td data-bbox="510 488 880 878">Over-Write Original Sample Files checkbox</td> <td data-bbox="880 488 1239 878"> replace the old files with the newly generated files. If you deselect this checkbox, a number is appended to the name for each newly generated file, and the original files are preserved. In BioLIMS mode, the file name must be unique. If the file name is not unique, then the sample file will not be written to the database. </td> </tr> </tbody> </table>	Choose...	To...	All Used Lanes radio button	generate a new Sample file for every lane marked Used.	Lanes marked for Extraction radio button (white markers)	generate a new Sample file only for each lane with a white lane marker.	Over-Write Original Sample Files checkbox	replace the old files with the newly generated files. If you deselect this checkbox, a number is appended to the name for each newly generated file, and the original files are preserved. In BioLIMS mode, the file name must be unique. If the file name is not unique, then the sample file will not be written to the database.
	Choose...	To...							
	All Used Lanes radio button	generate a new Sample file for every lane marked Used.							
Lanes marked for Extraction radio button (white markers)	generate a new Sample file only for each lane with a white lane marker.								
Over-Write Original Sample Files checkbox	replace the old files with the newly generated files. If you deselect this checkbox, a number is appended to the name for each newly generated file, and the original files are preserved. In BioLIMS mode, the file name must be unique. If the file name is not unique, then the sample file will not be written to the database.								
All Used Lanes radio button	generate a new Sample file for every lane marked Used.								
Lanes marked for Extraction radio button (white markers)	generate a new Sample file only for each lane with a white lane marker.								
Over-Write Original Sample Files checkbox	replace the old files with the newly generated files. If you deselect this checkbox, a number is appended to the name for each newly generated file, and the original files are preserved. In BioLIMS mode, the file name must be unique. If the file name is not unique, then the sample file will not be written to the database.								

To extract the data without changing current tracker lines: *(continued)*

Step	Action																
<p data-bbox="400 183 417 204">3</p>	<p data-bbox="471 183 1103 240">In the Project File portion of the dialog box you can take the following action:</p>																
	<table border="1"> <thead> <tr> <th data-bbox="467 285 830 321">Choose...</th> <th data-bbox="830 285 1174 321">To...</th> </tr> </thead> <tbody> <tr> <td data-bbox="467 321 830 578"> <p data-bbox="481 329 783 386">Add Sample Files to Project checkbox</p> </td> <td data-bbox="830 321 1174 578"> <ul style="list-style-type: none"> <li data-bbox="848 329 1150 443">◆ Create a new project, choose Create a New Project from the pop-up menu. <li data-bbox="848 459 1150 573">◆ Use an open project, choose Use the Open Project from the pop-up menu. </td> </tr> <tr> <td data-bbox="467 578 830 834"> <p data-bbox="481 586 770 643">Auto-Analyze New Sample Files checkbox</p> </td> <td data-bbox="830 578 1174 834"> <p data-bbox="848 586 1170 667">automatically analyze the new Sample files after extraction is finished.</p> <p data-bbox="848 691 1170 829">Deselect this if you want the Sample files extracted but not analyzed. This option is only available if the Sample files are being added to a project.</p> </td> </tr> <tr> <td data-bbox="467 834 830 902"> <p data-bbox="481 846 796 870">Analyze All Files radio button</p> </td> <td data-bbox="830 834 1174 902"> <p data-bbox="848 846 1143 902">analyze all the new Sample files created from the gel.</p> </td> </tr> <tr> <td data-bbox="467 902 830 971"> <p data-bbox="481 914 729 938">Print Results checkbox</p> </td> <td data-bbox="830 902 1174 971"> <p data-bbox="848 914 1130 971">print the results for all new Sample files after analysis.</p> </td> </tr> <tr> <td data-bbox="467 971 830 1065"> <p data-bbox="481 982 783 1039">Use Sample Sheet Settings checkbox</p> </td> <td data-bbox="830 971 1174 1065"> <p data-bbox="848 982 1157 1065">analyze and print the Sample files as designated in the Sample Sheet.</p> </td> </tr> <tr> <td data-bbox="467 1065 830 1333"> <p data-bbox="481 1076 756 1133">Save Gel After Extraction checkbox</p> </td> <td data-bbox="830 1065 1174 1333"> <p data-bbox="848 1076 1170 1157">save tracker lines and other gel file settings to the gel file after extracting Sample file data.</p> <p data-bbox="848 1182 1170 1320">If you do not select this option, the setting used for the extraction are discarded when you close the gel file without saving.</p> </td> </tr> <tr> <td data-bbox="360 1341 458 1409"> <p data-bbox="400 1349 417 1370">4</p> </td> <td data-bbox="458 1341 1184 1409"> <p data-bbox="471 1349 1116 1406">When all of the information in the Extract Lanes dialog box is correct, click OK to begin extracting data.</p> </td> </tr> </tbody> </table>	Choose...	To...	<p data-bbox="481 329 783 386">Add Sample Files to Project checkbox</p>	<ul style="list-style-type: none"> <li data-bbox="848 329 1150 443">◆ Create a new project, choose Create a New Project from the pop-up menu. <li data-bbox="848 459 1150 573">◆ Use an open project, choose Use the Open Project from the pop-up menu. 	<p data-bbox="481 586 770 643">Auto-Analyze New Sample Files checkbox</p>	<p data-bbox="848 586 1170 667">automatically analyze the new Sample files after extraction is finished.</p> <p data-bbox="848 691 1170 829">Deselect this if you want the Sample files extracted but not analyzed. This option is only available if the Sample files are being added to a project.</p>	<p data-bbox="481 846 796 870">Analyze All Files radio button</p>	<p data-bbox="848 846 1143 902">analyze all the new Sample files created from the gel.</p>	<p data-bbox="481 914 729 938">Print Results checkbox</p>	<p data-bbox="848 914 1130 971">print the results for all new Sample files after analysis.</p>	<p data-bbox="481 982 783 1039">Use Sample Sheet Settings checkbox</p>	<p data-bbox="848 982 1157 1065">analyze and print the Sample files as designated in the Sample Sheet.</p>	<p data-bbox="481 1076 756 1133">Save Gel After Extraction checkbox</p>	<p data-bbox="848 1076 1170 1157">save tracker lines and other gel file settings to the gel file after extracting Sample file data.</p> <p data-bbox="848 1182 1170 1320">If you do not select this option, the setting used for the extraction are discarded when you close the gel file without saving.</p>	<p data-bbox="400 1349 417 1370">4</p>	<p data-bbox="471 1349 1116 1406">When all of the information in the Extract Lanes dialog box is correct, click OK to begin extracting data.</p>
	Choose...	To...															
	<p data-bbox="481 329 783 386">Add Sample Files to Project checkbox</p>	<ul style="list-style-type: none"> <li data-bbox="848 329 1150 443">◆ Create a new project, choose Create a New Project from the pop-up menu. <li data-bbox="848 459 1150 573">◆ Use an open project, choose Use the Open Project from the pop-up menu. 															
	<p data-bbox="481 586 770 643">Auto-Analyze New Sample Files checkbox</p>	<p data-bbox="848 586 1170 667">automatically analyze the new Sample files after extraction is finished.</p> <p data-bbox="848 691 1170 829">Deselect this if you want the Sample files extracted but not analyzed. This option is only available if the Sample files are being added to a project.</p>															
	<p data-bbox="481 846 796 870">Analyze All Files radio button</p>	<p data-bbox="848 846 1143 902">analyze all the new Sample files created from the gel.</p>															
	<p data-bbox="481 914 729 938">Print Results checkbox</p>	<p data-bbox="848 914 1130 971">print the results for all new Sample files after analysis.</p>															
<p data-bbox="481 982 783 1039">Use Sample Sheet Settings checkbox</p>	<p data-bbox="848 982 1157 1065">analyze and print the Sample files as designated in the Sample Sheet.</p>																
<p data-bbox="481 1076 756 1133">Save Gel After Extraction checkbox</p>	<p data-bbox="848 1076 1170 1157">save tracker lines and other gel file settings to the gel file after extracting Sample file data.</p> <p data-bbox="848 1182 1170 1320">If you do not select this option, the setting used for the extraction are discarded when you close the gel file without saving.</p>																
<p data-bbox="400 1349 417 1370">4</p>	<p data-bbox="471 1349 1116 1406">When all of the information in the Extract Lanes dialog box is correct, click OK to begin extracting data.</p>																

How the GeneScan Analysis Software Names Sample Files

Naming Process The following table lists the process the GeneScan Analysis Software follows when naming generated Sample files:

If the generated Sample file...	Then GeneScan uses the...
has an associated file name in the Sample Sheet	file name in the Sample Sheet.
does not have an associated file name, but has a sample name ^a	lane number of the sample concatenated in front of the sample name, for example 02_Test1.
does not have an associated file name, or a sample name ^a	lane number of the sample concatenated in front of the name Sample file, for example 02_Sample File.
has the same name as a previously generated Sample file	lane number of the sample in parentheses concatenated to the file name, for example Test1.1, Test2.2.

a. These defaults can be edited in GS Collection. Refer to the *ABI PRISM 377 Instrument User's Manual*.

Saving Gel Files After Editing Tracking

Introduction The Gel File window keeps track of whether or not you have edited data, and when appropriate, will enable the Save command in the File menu. You may save all the information, or selected parts of the gel file.

Saving Files After Editing Tracking **IMPORTANT** If you save changes to the gel file, the original tracking information is overwritten. You can retrieve the originally calculated tracking only by choosing Track or Track & Extract Gel from the Gel menu to retrack the gel. See "Tracking Lanes and Extracting Data" on page 2-59.

You have the following options:

If you...	Then...
choose the Save command	The contents of the Gel File window are saved to the gel file.
select the checkbox labeled Save Gel after Extraction in the Extract Lane dialog box	It is not necessary to use the Save command to save the gel file.

For information regarding saving selected portions of gel files and archiving gel files, see "Saving Gel Files" on page 8-4.

Analyzing Sample Files

3

Introduction

In This Chapter Topics in this chapter include the following:

Topics	See page
About Sample Files	3-2
Opening Sample Files	3-4
About The Sample File Window	3-5
Sample Results View	3-6
Sample Info View	3-8
Size Curve View	3-13
Raw Data View	3-15
EPT Data View	3-17
Analyzing a Sample File	3-19

About Sample Files

What Sample Files Contain

Sample files contain data generated from the following:

Generated from...	Using...						
lane information from gel files	ABI™ 373 or ABI PRISM® 377 instruments. After extracting lanes from a gel file, Sample file data can be analyzed as follows:						
	<table border="1"> <thead> <tr> <th>Analyze Sample file data...</th> <th>For information, see...</th> </tr> </thead> <tbody> <tr> <td>directly from the Sample file</td> <td>“Analyzing a Sample File” on page 3-19.</td> </tr> <tr> <td>from the Analysis Control window in a project</td> <td>“Analyzing Sample Files: Using the Analysis Control Window” on page 5-6.</td> </tr> </tbody> </table>	Analyze Sample file data...	For information, see...	directly from the Sample file	“Analyzing a Sample File” on page 3-19.	from the Analysis Control window in a project	“Analyzing Sample Files: Using the Analysis Control Window” on page 5-6.
	Analyze Sample file data...	For information, see...					
directly from the Sample file	“Analyzing a Sample File” on page 3-19.						
from the Analysis Control window in a project	“Analyzing Sample Files: Using the Analysis Control Window” on page 5-6.						
capillary electrophoresis	ABI PRISM® 310 instrument.						

How GeneScan Generates Sample Files

The GeneScan® Analysis Software creates Sample files after extracting lanes. The information from each lane in the gel file is tracked and extracted, and the resulting Sample files are placed in their respective sample folder. If you change tracking, lane assignment, or Sample Sheet information, you have to regenerate the Sample files.

The software consults Sample Sheet information to determine whether a lane is used (contains sample). The lane tracker uses this information to assign lane numbers to the tracker lines. In addition, the GeneScan Analysis Software only extracts those lanes identified as Used.

Ways to Generate Sample Files

There are two ways to generate sample files:

Software	Instrument	Generates
Data Collection software	ABI PRISM 310	Sample file for each injection.
GeneScan Analysis Software	ABI 373 or ABI PRISM 377	Sample file for each lane.

How GeneScan Analyzes Sample Files

The GeneScan Analysis Software performs the following steps in analyzing Sample files:

Step	Action
1	Processes the raw data signals to generate analyzed data signal and then uses the analyzed signals to detect the signal peaks associated with DNA fragments.
2	Performs size calling by identifying the peaks of the in-lane size standard found in each sample.
3	<p>Determines the fragment size of each experimental peak within the sample based on the size calling curve generated using the size standard peaks, the selected size calling method, and by comparing it to the pre-defined size standard file.</p> <p>The algorithmic steps of the process from raw data to analyzed data are as follows:</p> <ul style="list-style-type: none">◆ Multicomponenting.◆ Baselineing.◆ Smoothing, if any.◆ Peak detection.

Opening Sample Files

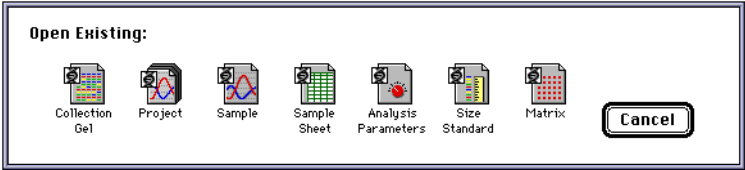
Introduction Sample files can be opened as separate files outside of projects, and display related information about each Sample file.

If you are interested in...	Then...
one or two Sample files	it is often more convenient to open Sample files individually and analyze or view the data without or opening an entire project.
multiple Sample files	use a project.

For information on opening Sample files from within:

- ◆ Projects, see “Accessing Sample Files” on page 5-6.
- ◆ BioLIMS™, see “Switching Between Sample File and BioLIMS Mode” on page E-10.

Procedure To open a Sample file as a separate file:

Step	Action
1	Choose Open from the File menu. The Open Existing dialog box appears. Note You can also double-click the Sample file name in the Finder. If the GeneScan Analysis Software is not running, the software starts and opens the Sample file.
	
2	Click the Sample icon. A directory dialog box appears.
3	In the dialog box, find and select the Sample file that you want to open.
4	Click Open. The Sample File window appears. See “About The Sample File Window” on page 3-5.

About The Sample File Window

What it Displays You can use the different display modes in the Sample File window to review the analyzed and raw data, and all pertinent data collection, sizing and Sample description information from a single window. The Sample Results view appears as the default.


Five Views The five views of the Sample File window are:

View	See page
Sample Results view	3-6
Sample Info view	3-8
Size Curve view	3-13
Raw Data view	3-15
EPT Data view	3-17

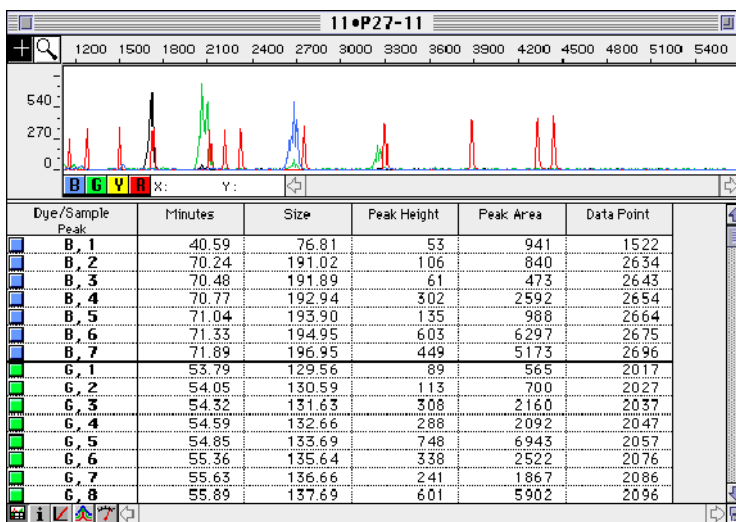
Sample Results View

What it Displays The Sample Results View displays the Sample file's analyzed data in both electropherogram and tabular data form.

Displaying the View The following table lists ways to display the Sample Results View:

To display the view...	Do this...
from the Sample file window	<ul style="list-style-type: none"> ◆ Click the button for the Sample Results view at the bottom left of the Sample file window,  or, ◆ Choose Sample Results (⌘ E) from the Sample menu.
from a project window	<ol style="list-style-type: none"> a. Select a sample or multiple samples in the Analysis Control or the Results Control window. b. Choose Sample Results from the Sample menu.

Sample Results View Example The following is an example of the Sample Results view:



Description of Columns

The following table describes the columns in the above figure:

This column	Identifies
Dye/Sample Peak	◆ Dye color. ◆ Peak number.
Minutes	The time, in minutes, from the start of the run to the time the fragment was detected.
Size	The number of base pairs in the fragment. This value is calculated automatically only if you: ◆ Run the size standard in the same lane or injection as the sample, and ◆ Perform size calling.
Peak Height	Signal size.
Peak Area	Area of the detected peak.
Data point	Data point of the fragment at its maximum peak height. Scan number for ABI 373, ABI 373XL, ABI PRISM 377 and ABI PRISM 377XL data.

Differences From the Results Display

The Sample Results view displays the same electropherogram and tabular data as the Results Display, with the following differences:


- ◆ One Sample file displayed.
 - ◆ Show or hide dye/sample data by clicking the buttons below the electropherogram.
 - ◆ Cannot display legends.
 - ◆ Cannot use custom plot colors.
-
-

Sample Info View

What it Displays Displays the following Sample file information:

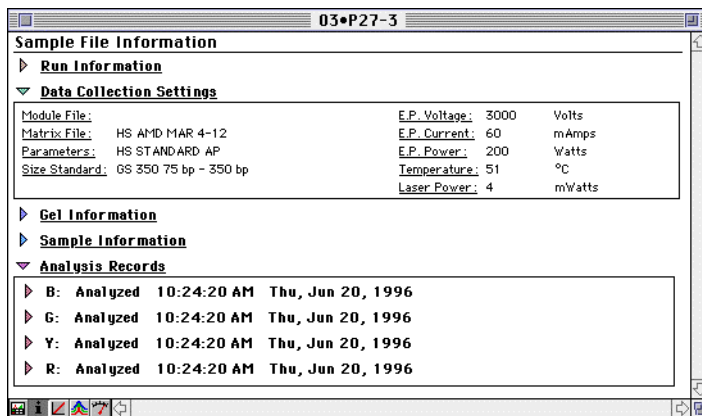
- ◆ Run and data collection information.
 - ◆ Gel information.
 - ◆ Sample information—this information can be edited.
 - ◆ Analysis records.
-

Displaying the View The following table lists ways to display the Sample Info View:

To display the view...	Do this...
from the Sample file window	<ul style="list-style-type: none">◆ Click the button for the Sample Info view at the bottom left of the Sample file window,  or,◆ Choose Sample Info (⌘ I) from the Sample menu.
from a project window	<ol style="list-style-type: none">a. Select a sample or multiple samples in the Analysis Control or the Results Control window.b. Choose Sample Info from the Sample menu. <p>The information is organized in five panels.</p> <p>Click the triangles to expand or collapse the panels to display specific information.</p>

Sample Info View Example

The following is an example of the Sample File window in Sample Info View:



Description of Information

The following tables list the information in the Sample Info View:

Run Information

Information found under the header	Information inserted from
<ul style="list-style-type: none"> ◆ User name. ◆ Instrument. ◆ Data Collection software version. ◆ Run date and start time. ◆ Run duration. ◆ Total data points collected. 	<ul style="list-style-type: none"> ◆ ABI PRISM 310 data-collection Run file and run information. ◆ ABI 373 or ABI PRISM 377 gel file (data collection Run file and run information). <p>This information is embedded in the gel file.</p>

Data Collection Settings

Information found under the header	Information is inserted from
<ul style="list-style-type: none"> ◆ Module File. ◆ Matrix File. ◆ Analysis Parameters. ◆ Size Standard. ◆ Electrophoresis voltage, current, and power (ABI 373 and ABI PRISM 377 only). ◆ Run Voltage, Injection Voltage, and Injection Duration (ABI PRISM 310 only). ◆ Temperature. ◆ Laser power. 	<p>ABI 373 or ABI PRISM 377 gel file (data collection Run File and run information).</p> <p>This information is embedded in the gel file.</p> <p>ABI PRISM 310 data collection Run file and run information.</p>

Gel Information (Polymer)

Information found under the header	Information is inserted from
ABI PRISM 310	
◆ Gel type.	Optional (user entered).
◆ Length to detector.	Data Collection information.
◆ Lot # and expiration date.	
ABI 373 and ABI PRISM 377	
◆ Gel type.	Optional (user entered).
◆ Name of gel file.	<p>Gel file (data collection Run file).</p> <p>This information is embedded in the gel file.</p>
◆ Gel percentage and thickness.	Optional (user entered).
◆ Well-to-read (separation) distance.	
◆ Number of channels.	Gel file (data collection Run file).
◆ Number of Lanes.	This information is embedded in the gel file.

Information found under the header	Information is inserted from										
<p>◆ Channel Averaging.</p> <p>Note Zero (0) indicates use of pre-averaging offscale data.</p> <p>The following is an example of how pre-averaging offscale data appears:</p> <table border="1" data-bbox="417 386 821 768"> <thead> <tr> <th data-bbox="424 391 669 427">Extraction method</th> <th data-bbox="669 391 814 427">Displayed</th> </tr> </thead> <tbody> <tr> <td data-bbox="424 427 669 467">3 channel averaging</td> <td data-bbox="669 427 814 467">3</td> </tr> <tr> <td data-bbox="424 467 669 565">3 channel averaging with pre-averaging offscale</td> <td data-bbox="669 467 814 565">30</td> </tr> <tr> <td data-bbox="424 565 669 634">3 channel averaging-weighted</td> <td data-bbox="669 565 814 634">- 3</td> </tr> <tr> <td data-bbox="424 634 669 768">3 channel averaging-weighted, pre-averaging offscale</td> <td data-bbox="669 634 814 768">- 30</td> </tr> </tbody> </table>	Extraction method	Displayed	3 channel averaging	3	3 channel averaging with pre-averaging offscale	30	3 channel averaging-weighted	- 3	3 channel averaging-weighted, pre-averaging offscale	- 30	Gel processing parameters.
Extraction method	Displayed										
3 channel averaging	3										
3 channel averaging with pre-averaging offscale	30										
3 channel averaging-weighted	- 3										
3 channel averaging-weighted, pre-averaging offscale	- 30										
◆ Range of data points extracted.											

Sample Information

Information found under header	Information is inserted from
Sample data and comment for each dye color.	<p>◆ ABI PRISM 310 Sample Sheet.</p> <p>◆ ABI 373 or ABI PRISM 377 gel file (Sample Sheet).</p> <p>This information is embedded in the gel file.</p>

Analysis Records

Information found under header	Information is inserted from
Date and time each color was analyzed, and more panel-display arrows.	Analysis information.


Dyes Within Analysis Records

Information found under header	Information is inserted from
◆ Analysis parameters file and range analyzed.	Analysis Settings.
◆ Whether baselined or multicomponented.	
◆ Data smoothing.	
◆ Peak detection threshold and minimum half-width.	
◆ Size standard file.	Analysis Control window.
◆ The dye color used for the standard, Sizing method, and range.	
◆ Standard peak detection threshold.	Analysis Settings.
◆ Split Peak correction.	
◆ Total number of peaks: found in sample & dye standard, defined in standard matched with standard peaks.	Analysis results.

Size Curve View

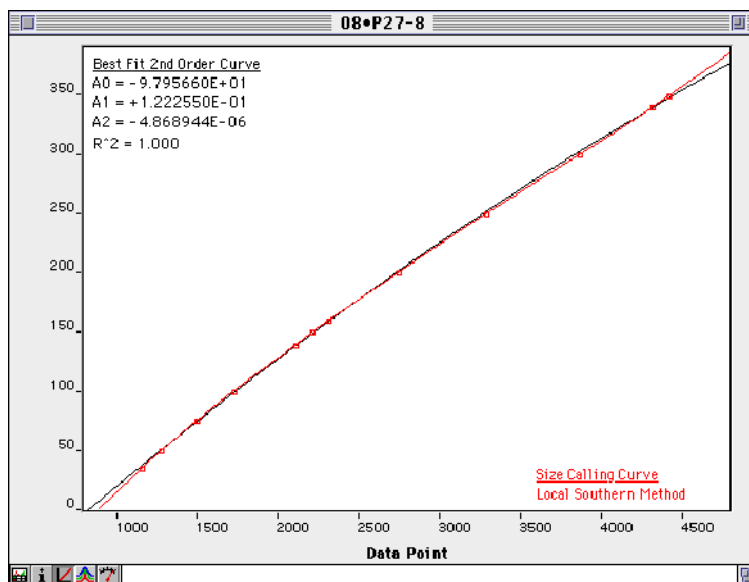
What it Displays Displays sizing curves for Sample files. The size curve is a measure of how well the internal size standard matches the standard definition, and whether or not it is linear.

Displaying the View The following table lists ways to display the Size Curve View:

To display the view...	Do this...								
from the Sample file window	<ul style="list-style-type: none"> ◆ Click the Size Curve button on the lower left of the Sample file window  or ◆ Select Size Curve (⌘ U) from the Sample menu. 								
from a project window	<p>a. Use the following table to select a sample or multiple samples in the Analysis Control or the Results Control window.</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th style="width: 50%;">If you want to...</th> <th style="width: 50%;">Then...</th> </tr> </thead> <tbody> <tr> <td>select a sample</td> <td>click the sample.</td> </tr> <tr> <td>you want to select several consecutive samples</td> <td>shift-click the first and last sample in the group you want to select.</td> </tr> <tr> <td>select several samples that are not next to each other</td> <td>⌘-click the samples.</td> </tr> </tbody> </table> <p>b. Choose Size Curve from the Sample menu. For each selected sample, a Sample file window opens and displays its Size Curve view.</p>	If you want to...	Then...	select a sample	click the sample.	you want to select several consecutive samples	shift-click the first and last sample in the group you want to select.	select several samples that are not next to each other	⌘-click the samples.
If you want to...	Then...								
select a sample	click the sample.								
you want to select several consecutive samples	shift-click the first and last sample in the group you want to select.								
select several samples that are not next to each other	⌘-click the samples.								

What the Size Curve Displays

The Size Curve view displays two curves, as shown in the figure below. For a description of the curves, see “Curves Described” below.



Curves Described

The following table describes the curves in the above figure:


Note Sizing errors due to anomalous mobilities are displayed as non-linear.

This curve	Represents
Red curve	The size calling curve, based on the size calling method used to analyze the data.
Black curve	<p>The best-fit least squares curve, which the GeneScan Analysis Software calculates for all samples, regardless of the size calling method.</p> <p>This curve is provided to help evaluate the linearity of the sizing curve.</p> <p>When the sizing curve and best-fit curve match, they overlap so you see only the size curve.</p>

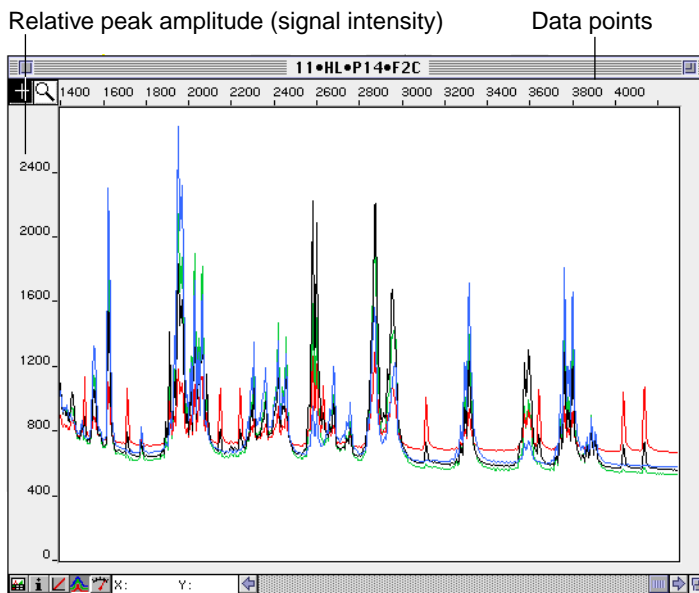
Raw Data View

What it Displays Displays the raw data collected for a sample. Raw data is non-baselined and non-multicomponented. This information is stored in the Sample file. The gel file does not need to be available.

Displaying the View The following table lists ways to display the Raw Data View:

To display the view...	Do this...
from the Sample file window	<ul style="list-style-type: none">◆ Click the button for the Raw Data view at the bottom left of the Sample file window,  or,◆ Choose Raw Data (⌘ R) from the Sample menu.
from a project window	<ol style="list-style-type: none">a. Select a sample or multiple samples in the Analysis Control or the Results Control window.b. Choose Raw Data (⌘ R) from the Sample menu.

Raw Data Example The following is an example of the Sample File window in Raw Data View:



What to Evaluate Use the Raw Data view to evaluate:


- ◆ Problems or noise in the baseline that could result in poor size calling.
- ◆ Areas with a lower signal may indicate bad tracking.
- ◆ Start and stop points for analysis.

For information on changing the horizontal (page 7-37) or vertical scale of the data (see page 7-38).

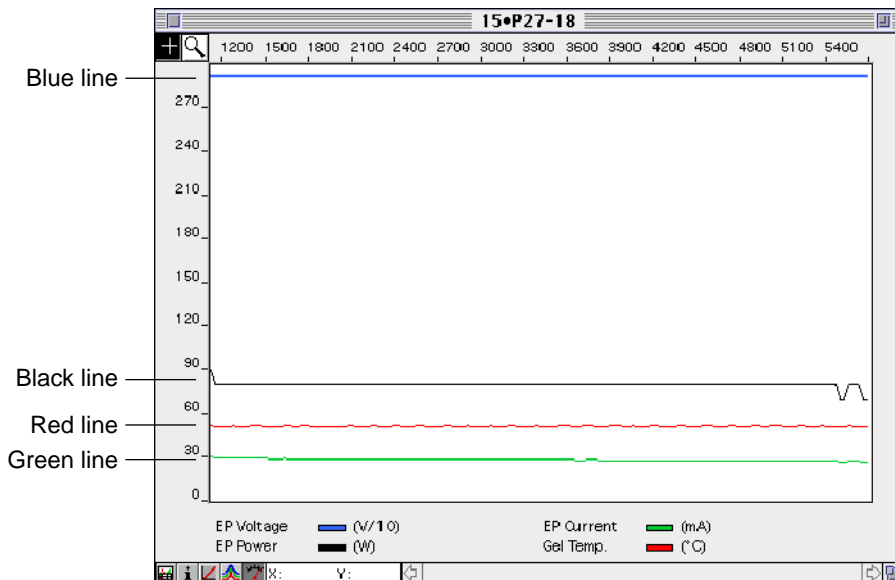
EPT Data View

What it Displays Displays the electrophoresis power and temperature information collected for a sample. This information is stored in the Sample file. The gel file does not need to be available.

Displaying the View The following table lists ways to display the EPT Data View:

To display the view....	Do this...
from a Sample file window	<ul style="list-style-type: none">◆ Click the button for the EPT Data view at the bottom left of the Sample file window,  or◆ Choose EPT Data (⌘ M) from the Sample menu.
from a project window	<ul style="list-style-type: none">a. Select a sample or multiple samples in the Analysis Control or the Results Control window.b. Select EPT Data from the Sample menu, or click the EPT data button.

EPT Data Example The following is an example of the Sample File window in EPT Data View:



Colored Lines Described The following table describes the lines in the above figure:

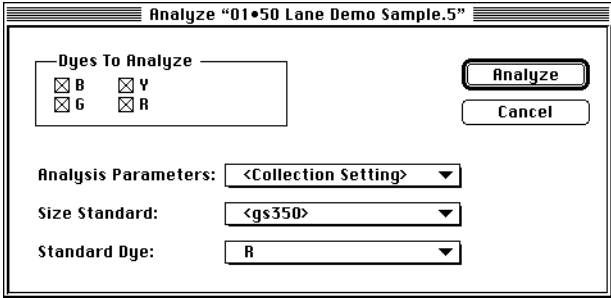
Line	Description
Blue	Electric voltage in volts/10
Black	Electric power in watts
Red	Gel temperature in °C
Green	Electric current in mA (mamp)

Analyzing a Sample File

Introduction The GeneScan Analysis Software analyzes raw data stored in Sample files according to parameters and standards that you select. You can use the analyzed data to detect peaks associated with DNA fragments, and identify those peaks with an established size standard.

Procedure

To analyze a sample file:

Step	Action										
1	<p>Choose Analyze "Sample File Name" from the Sample menu (⌘ Y). The Analyze Sample file dialog box appears.</p> 										
2	<p>From the Analyze Sample file dialog box, select one of the following options:</p> <table border="1"> <thead> <tr> <th>Choose...</th> <th>To select...</th> </tr> </thead> <tbody> <tr> <td>Dyes to Analyze checkboxes</td> <td>any number of dyes to analyze.</td> </tr> <tr> <td>Analyze Parameters pop-up menu</td> <td>from the default parameters or any parameter files in the folder location specified in the application preferences.</td> </tr> <tr> <td>Size Standard pop-up menu</td> <td>from the default standard, or any standard files in the folder location that you specify in the application preferences.</td> </tr> <tr> <td>Standard Dye pop-up menu</td> <td>the inlane standard dye.</td> </tr> </tbody> </table>	Choose...	To select...	Dyes to Analyze checkboxes	any number of dyes to analyze.	Analyze Parameters pop-up menu	from the default parameters or any parameter files in the folder location specified in the application preferences.	Size Standard pop-up menu	from the default standard, or any standard files in the folder location that you specify in the application preferences.	Standard Dye pop-up menu	the inlane standard dye.
Choose...	To select...										
Dyes to Analyze checkboxes	any number of dyes to analyze.										
Analyze Parameters pop-up menu	from the default parameters or any parameter files in the folder location specified in the application preferences.										
Size Standard pop-up menu	from the default standard, or any standard files in the folder location that you specify in the application preferences.										
Standard Dye pop-up menu	the inlane standard dye.										

To analyze a sample file: *(continued)*

Step	Action
3	After analysis, evaluate the results. For information on evaluating the results, see Chapter 7, "Evaluating Analysis Results."

Installing a New Matrix File

Use the following procedure to install a new matrix file for the Sample file that you want to analyze.

For information on attaching the new matrix to an ABI 373 or ABI PRISM 377 gel file, see "Installing New Matrix Information" on page 2-24.

To install a new matrix file:

Step	Action
1	Choose Install New Matrix from the Sample menu. A directory dialog box appears. The Folder Preferences settings determine where the GeneScan Analysis Software looks for the matrix file. For more information, see "Defining Folder Locations" on page 5-41.
2	Select the new matrix file in the dialog box and click Open. A message appears when the matrix is successfully assigned.
3	Re-Analyze the Sample file. Applying a new matrix file clears previous analysis information, so you must re-analyze the file.

4

Creating a Project

Introduction

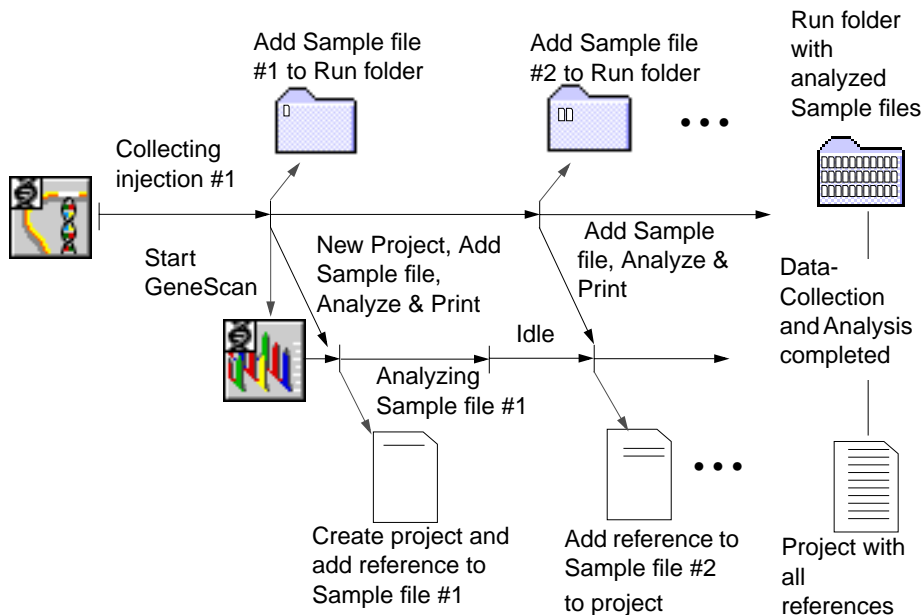
In This Chapter Topics in this chapter include the following:

Topics	See page
Automatic Analysis and Project Creation Process	4-2
Setting Up for Automatic Analysis	4-4
Using a Project to Manage Sample Files	4-8
Working with Project Files	4-10
Finding Missing Sample Files	4-14

Automatic Analysis and Project Creation Process

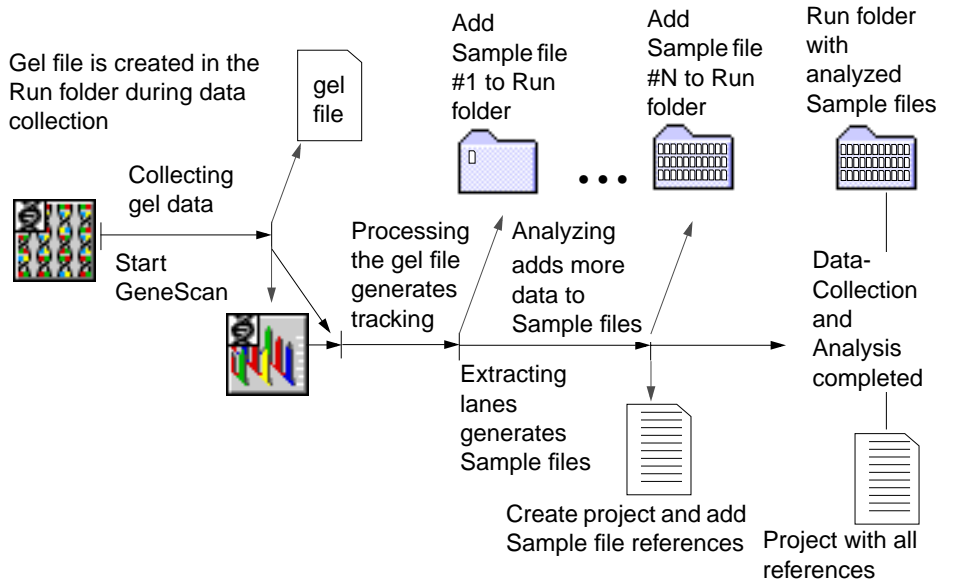
Introduction The following two diagrams illustrate the process of automatic analysis and project creation for data from the ABI PRISM® 310 and from the ABI™ 373 and ABI PRISM® 377 instruments.

Process Using the ABI PRISM 310 The following diagram illustrates data analysis using the ABI PRISM 310 instrument.



**Process Using the
ABI 373 and the
ABI PRISM 377**

The following diagram illustrates data analysis using the ABI 373 and the ABI PRISM 377 instruments.



Setting Up for Automatic Analysis

Introduction To set up the GeneScan® Analysis Software for automatic analysis after data-collection, you must have previously defined analysis parameters and size standards.

For more information see:

- ◆ “Defining Analysis Parameters” on page 5-18.
- ◆ “Defining the Size Standard” on page 5-31.

Procedure

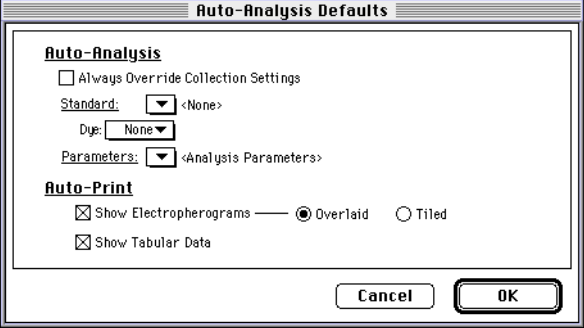
To set up for automatic analysis after data collection:

Step	Action
Complete the following steps in the data collection software.	
1	Set the GeneScan Run default preferences to auto-analyze and use the pop-up menu to locate and select the GeneScan Analysis Software.
2	<p>In the GeneScan Sample Sheet, for each sample to be analyzed do the following:</p> <ul style="list-style-type: none">◆ Enter the sample name. This field must be completed for the samples to be active in the Injection or Run Sheet.◆ Indicate which dye is the standard.◆ Select the checkbox labeled Pres (Present) for each dye/sample you want auto-analyzed.◆ Select any additional checkboxes. <p>IMPORTANT If you plan to use the Genotyper® software, then you must complete the Sample Info box correctly. For more information, refer to the <i>Genotyper User's Manual</i>.</p>

To set up for automatic analysis after data collection: *(continued)*

Step	Action						
3	<p>In the Injection List or Run Sheet for each applicable sample, select AutoAnalyze, and choose from the appropriate pop-up menus one of the following:</p> <ul style="list-style-type: none"> ◆ Matrix file. ◆ Analysis parameters. ◆ Size standard. <p>The following table lists considerations when choosing a size standard.</p> <table border="1" data-bbox="577 505 1239 919"> <thead> <tr> <th data-bbox="577 505 907 544">If you choose a...</th> <th data-bbox="907 505 1239 544">Then GeneScan...</th> </tr> </thead> <tbody> <tr> <td data-bbox="577 544 907 732">Size Standard definition file in the Collection Injection List or Run Sheet, but you do not specify which dye is the standard in the Sample Sheet</td> <td data-bbox="907 544 1239 732">performs analysis using the dye specified in the Auto-Analysis defaults.</td> </tr> <tr> <td data-bbox="577 732 907 919">dye as the standard in the Sample Sheet, but do not specify a size standard definition file in the Collection Injection List or Run Sheet</td> <td data-bbox="907 732 1239 919">does not perform size calling.</td> </tr> </tbody> </table>	If you choose a...	Then GeneScan...	Size Standard definition file in the Collection Injection List or Run Sheet, but you do not specify which dye is the standard in the Sample Sheet	performs analysis using the dye specified in the Auto-Analysis defaults.	dye as the standard in the Sample Sheet, but do not specify a size standard definition file in the Collection Injection List or Run Sheet	does not perform size calling.
If you choose a...	Then GeneScan...						
Size Standard definition file in the Collection Injection List or Run Sheet, but you do not specify which dye is the standard in the Sample Sheet	performs analysis using the dye specified in the Auto-Analysis defaults.						
dye as the standard in the Sample Sheet, but do not specify a size standard definition file in the Collection Injection List or Run Sheet	does not perform size calling.						
4	Select Auto Print to print automatically.						
Complete the following steps in the GeneScan Analysis Software:							
1	<p>Take the following action for the ABI 373 and the ABI PRISM 377 data:</p> <ol style="list-style-type: none"> a. Start the GeneScan Analysis Software. b. From the Gel menu and the Auto-Processing submenu, ensure that Auto-Track Gel and Extract Lanes after Auto-Tracking are selected. <p>IMPORTANT Before choosing Auto-Track Lane, verify that the comb type is set correctly by choosing Preferences from the Settings menu and Gel Preferences from the submenu.</p> <p>For more information, see “How to Set Gel Processing Preferences” on page 2-3.</p>						

To set up for automatic analysis after data collection: *(continued)*

Step	Action
<p>2</p>	<p>If you are using the Auto-Analysis defaults, choose Auto-Analysis Defaults from the Settings menu.</p> <p>The Auto-Analysis Defaults dialog box appears.</p>  <p>The parameters set in the Auto-Analysis Defaults dialog box apply when you have:</p> <ul style="list-style-type: none"> ◆ Specified <Analysis Defaults> in the data-collection software. ◆ Selected the checkbox labeled Always Override Collection Settings.
<p>3</p>	<p>Select the checkbox labeled Always Override Collection Settings to have the parameters set take precedence over the following files specified in the Data Collection software:</p> <ul style="list-style-type: none"> ◆ Size standard. ◆ Analysis parameters. ◆ Dye standard.
<p>4</p>	<p>Choose a new Size Standard if:</p> <ul style="list-style-type: none"> ◆ <Analysis Defaults> is specified for the size standard in the data collection settings, or ◆ To override the specified size standard.
<p>5</p>	<p>Choose the Dye pop-up menu to specify the dye that represents the internal size standard you are running with the samples.</p>
<p>6</p>	<p>Choose the analysis parameters you want to use from the Parameters pop-up menu.</p> <p>To use the default parameters, choose <Analysis Parameters>.</p>

To set up for automatic analysis after data collection: *(continued)*

Step	Action						
7	In the Auto-Print section choose, you have the following options:						
	<table border="1"> <thead> <tr> <th data-bbox="510 251 873 284">Choose...</th> <th data-bbox="873 251 1229 284">To print...</th> </tr> </thead> <tbody> <tr> <td data-bbox="510 284 873 576">Show Electropherograms</td> <td data-bbox="873 284 1229 576"> electropherograms. Select the appropriate radio button to specify whether the electropherograms for the four dyes appear: ♦ Together in one panel (overlaid) or in ♦ Separate panels (tiled). </td> </tr> <tr> <td data-bbox="510 576 873 625">Show Tabular data</td> <td data-bbox="873 576 1229 625">tabular data.</td> </tr> </tbody> </table>	Choose...	To print...	Show Electropherograms	electropherograms. Select the appropriate radio button to specify whether the electropherograms for the four dyes appear: ♦ Together in one panel (overlaid) or in ♦ Separate panels (tiled).	Show Tabular data	tabular data.
	Choose...	To print...					
Show Electropherograms	electropherograms. Select the appropriate radio button to specify whether the electropherograms for the four dyes appear: ♦ Together in one panel (overlaid) or in ♦ Separate panels (tiled).						
Show Tabular data	tabular data.						
8	<p>Click OK.</p> <p>The GeneScan Analysis Software is now prepared to perform automatic analysis.</p> <p>Note The settings that you specify are initial settings. Make two or three trial runs, fine tuning the parameters with each run, to determine which parameters work the best for a particular protocol.</p>						

Using a Project to Manage Sample Files

What is a Project A project is a file containing references to a set of Sample files that you want to analyze and display together. The project contains Analysis Control and Results Control windows that allow you to analyze specific dye/samples and display the results of analysis.

Why Create a Project You can create a new project and add any combination of Sample files, allowing you to analyze and display samples from different runs.

Adding a Sample file to the project sets up a link between the project and the Sample file. The file itself is not imported into the project.

For more information, see “Creating a New Project” on page 4-10.

When GeneScan Creates Projects The GeneScan Analysis Software creates a project file when Sample files are created during automatic analysis.

Note When the GeneScan Analysis Software automatically creates a project, the default file name is the same as the gel file.

The following table lists when the GeneScan Analysis Software creates a project file:

When using this instrument...	A project is created...
ABI PRISM 310	immediately after the first sample run. Note The default project name is GeneScan Project - “date”.
ABI 373 and ABI PRISM 377	when the software extracts the lanes of the gel file to create the Sample files. Note The default project name is the same as the gel with the following extension added to the end, “.π”.

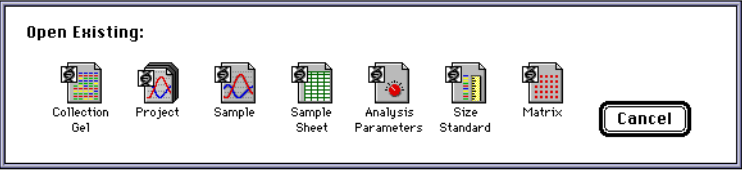
Where to Store Projects Keep the Sample files in the same location on the hard disk relative to the project file so the GeneScan Analysis Software can locate them when the project is opened.

If Sample files are moved and the program does not find them when you open the project, see “Finding Missing Sample Files” on page 4-14.

Working with Project Files

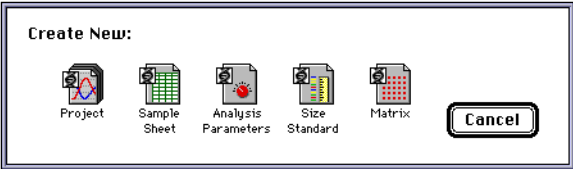
Opening an Existing Project

To open an existing project:

Step	Action
1	<p>Choose Open from the File menu.</p> <p>The Open Existing dialog box appears.</p> <p>For information on opening a Sample file from the BioLIMS™ database, see “Switching Between Sample File and BioLIMS Mode” on page E-10.</p>
	
2	<p>Click the Project icon.</p> <p>A directory dialog box appears.</p>
3	<p>In the dialog box, find and select the project you want to open.</p>
4	<p>Click Open.</p> <p>The project opens in an Analysis Control window.</p>

Creating a New Project

To create a new project:

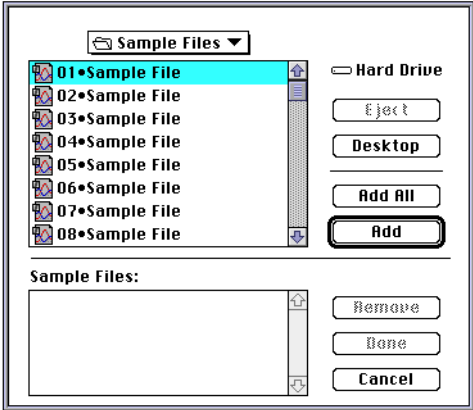
Step	Action
1	<p>Choose New from the File menu.</p> <p>The Create New dialog box appears.</p>
	

To create a new project: *(continued)*

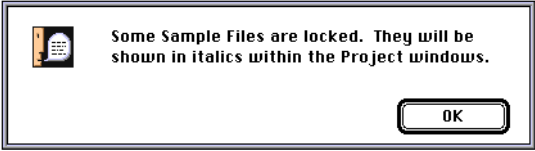
Step	Action						
2	<p>Click the Project icon.</p> <p>An untitled Analysis Control window appears.</p> <p>For information on using the Analysis Control window, see page 5-4.</p>						
3	<p>There are two ways to add files to a project:</p> <table border="1"> <thead> <tr> <th>To add...</th> <th>Choose...</th> </tr> </thead> <tbody> <tr> <td>sample files you select to the open project</td> <td> <p>Add Sample Files (⌘ B).</p> <p>The Add Sample dialog box appears (see “Using the Add Sample Dialog Box” below).</p> </td> </tr> <tr> <td>a sample currently open to the open project</td> <td>Add “file name”.</td> </tr> </tbody> </table>	To add...	Choose...	sample files you select to the open project	<p>Add Sample Files (⌘ B).</p> <p>The Add Sample dialog box appears (see “Using the Add Sample Dialog Box” below).</p>	a sample currently open to the open project	Add “file name”.
To add...	Choose...						
sample files you select to the open project	<p>Add Sample Files (⌘ B).</p> <p>The Add Sample dialog box appears (see “Using the Add Sample Dialog Box” below).</p>						
a sample currently open to the open project	Add “file name”.						

Using the Add Sample Dialog Box

To use the Add Sample dialog box:

Step	Action
1	<p>Find the folder containing the samples that you want to add from the top scroll box.</p> 

To use the Add Sample dialog box: *(continued)*

Step	Action								
2	You can take the following action:								
	<table border="1"> <thead> <tr> <th>If...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>you want to select a single Sample file</td> <td>double-click the file or select the file and click Add.</td> </tr> <tr> <td>you want to select all the Sample files</td> <td>click Add All.</td> </tr> <tr> <td>you want to add a random selection of Sample files</td> <td>double-click each file name.</td> </tr> </tbody> </table>	If...	Then...	you want to select a single Sample file	double-click the file or select the file and click Add.	you want to select all the Sample files	click Add All.	you want to add a random selection of Sample files	double-click each file name.
	If...	Then...							
	you want to select a single Sample file	double-click the file or select the file and click Add.							
you want to select all the Sample files	click Add All.								
you want to add a random selection of Sample files	double-click each file name.								
3	Click Done when you have added all the Sample files.								
	<table border="1"> <thead> <tr> <th>If the...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>Sample files appear in the Analysis Control window</td> <td>see "Analyzing Sample Files: Using the Analysis Control Window" on page 5-6.</td> </tr> <tr> <td>Locked files alert appears (see below)</td> <td>go to "Unlocking Sample Files" below.</td> </tr> </tbody> </table> <div style="border: 1px solid black; padding: 10px; margin-top: 10px;">  </div>	If the...	Then...	Sample files appear in the Analysis Control window	see "Analyzing Sample Files: Using the Analysis Control Window" on page 5-6.	Locked files alert appears (see below)	go to "Unlocking Sample Files" below.		
If the...	Then...								
Sample files appear in the Analysis Control window	see "Analyzing Sample Files: Using the Analysis Control Window" on page 5-6.								
Locked files alert appears (see below)	go to "Unlocking Sample Files" below.								

Unlocking Sample Files

If the Sample files added to a project are locked, the GeneScan Analysis Software does not allow changes to them. You cannot analyze locked files. When you add locked files, an alert appears.

To unlock locked files:

Step	Action
1	Save and close the project.
2	In the Finder, select the applicable Sample files.
3	Choose Get Info from the File menu.
4	In the lower left-hand corner of each Sample Info window, deselect the checkbox labeled locked.


To unlock locked files:

Step	Action
5	<p>Close the window and reopen the project.</p> <p>Note If you did not close the project before unlocking the files, once the files are unlocked, close and open the project so the GeneScan Analysis Software recognizes that the Sample files are unlocked.</p>

Removing Samples from a Project

Note A removed Sample file is not deleted from the hard disk. The reference is removed from the project.

To remove Sample files from a project:

Step	Action
1	<p>Select the file or files in the Analysis Control window or the Results Control window that you want to remove.</p> <p>Note Shift-click to select multiple consecutive file or ⌘-click to select multiple files that are not consecutive.</p>
2	<p>Choose Remove Sample Files from the Project menu, or press the Delete key.</p> <p>A warning dialog box appears.</p> <div data-bbox="521 847 1055 982" style="border: 1px solid black; padding: 10px; margin: 10px 0;"><p>1 Sample Files are selected. Are you sure you want to remove them?</p><p>Cancel Remove</p></div>
3	<p>Click Remove.</p>

Finding Missing Sample Files

When are Files Considered Lost The project and related Sample files are usually located in the Run folder created by the Data Collection software. If the Sample files or the project are moved so they are no longer in the same relative position, the GeneScan Analysis Software might not be able to locate the Sample files when the project is opened. Usually, this only occurs if the Sample files are moved to another disk drive, another server on a network, or to another disk partition on the hard drive.

When an Alert Appears If the GeneScan Analysis Software does not locate the Sample files associated with a project, an alert box appears and the Sample file names appear dimmed when the project opens.

Searching for Missing Sample Files You can re-establish the links between the Sample files and the project by choosing Find Missing Sample Files from the Project menu and choosing one of the following options from the submenu.

Note Version 2.1 and later versions of the software allows you to choose how you want to search for missing Sample files.

Searching for missing sample files:

Choose...	If you...	Description
Fast Search	suspect that the GeneScan Analysis Software could not find the missing Sample files because they are located on an unmounted external storage device, or diskette.	Once the volume is mounted, or the diskette containing the files is inserted, Fast Search finds them immediately.
Search a Folder	know what folder contains the missing Sample file.	Specify a folder in which to search for missing Sample files. The GeneScan Analysis Software then immediately locates the Sample files and re-establishes links to the project.

Searching for missing sample files: *(continued)*

Choose...	If you...	Description
Exhaustive Search	do not know where any of the specified missing Sample files are located.	The GeneScan Analysis Software searches all mounted disk drives, and available servers. When the files are found, the software re-establishes links to the project.

**Re-Establishing
Links**

To re-establish the links with Sample files:

Step	Action
1	Click the dimmed file name to select it. When the missing file is selected the Find "file name" command becomes active. The name of the missing file appears inside the quotation marks.
2	Choose Find "file name". A file dialog box appears.
3	In the file dialog box, locate the proper folder and file.
4	Click Open or double-click the file name.

Analyzing Project Files

5

Introduction

In This Chapter Topics in this chapter include the following:

Topics	See page
Analyzing Project Files: About the Analysis Control Window	5-2
Analyzing Sample Files: Using the Analysis Control Window	5-6
About the Analysis Parameters	5-17
Using Analysis Parameter Files	5-24
About Size Standards	5-29
Defining the Size Standard	5-31
Using Size Standards	5-36
Defining Folder Locations	5-41

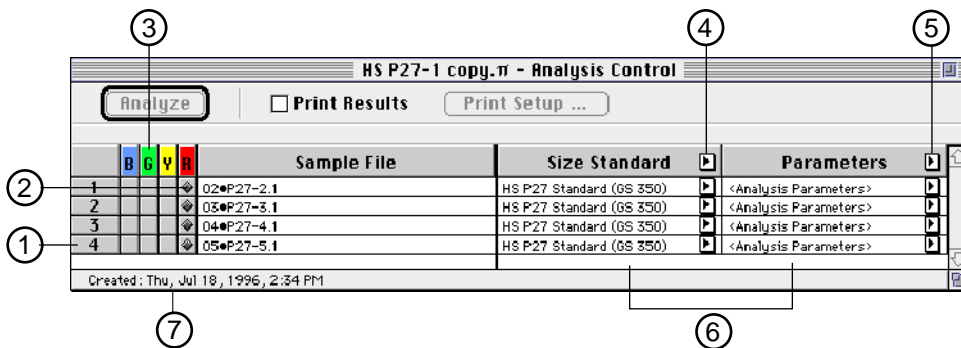
Analyzing Project Files: About the Analysis Control Window

Introduction When a project is opened the Analysis Control window appears. The Analysis Control window is the main window of a project. You can use this window to specify the following for each sample in the project:

- ◆ Dye that represents the size standard you ran with the sample.
- ◆ Size standard.
- ◆ Analysis parameters.
- ◆ Specific dyes to be analyzed.
- ◆ Optionally, format the document for printing and print the results automatically.

For more information, see “Analyzing Sample Files: Using the Analysis Control Window” on page 5-6.

Analysis Control Window Example The following is an example of the Analysis Control window:



**Analysis Control
Window
Description**

The following table describes the above diagram:

No.	Description						
1	Use these columns to choose the dye colors to analyze and to specify which is the size standard.						
2	Diamonds mark the standards. For more information, see "What are Size Standards" on page 5-29.						
3	⌘-click a dye/sample field to specify that dye/sample as the standard.						
4	<p>From this pop-up menu, you can choose:</p> <ul style="list-style-type: none"> ◆ <Collection Settings>. ◆ A user-defined size standard. ◆ Define a new size standard for that sample. <p>For more information, see:</p> <table border="1" data-bbox="514 688 1200 808"> <thead> <tr> <th data-bbox="514 688 1053 727">Topic</th> <th data-bbox="1053 688 1200 727">See page</th> </tr> </thead> <tbody> <tr> <td data-bbox="514 727 1053 766">About Size Standards</td> <td data-bbox="1053 727 1200 766">5-29</td> </tr> <tr> <td data-bbox="514 766 1053 808">Using Size Standards</td> <td data-bbox="1053 766 1200 808">5-36</td> </tr> </tbody> </table>	Topic	See page	About Size Standards	5-29	Using Size Standards	5-36
Topic	See page						
About Size Standards	5-29						
Using Size Standards	5-36						
5	<p>From this pop-up menu, choose an analysis parameters file.</p> <p>For more information, see:</p> <table border="1" data-bbox="514 938 1200 1058"> <thead> <tr> <th data-bbox="514 938 1053 977">Topic</th> <th data-bbox="1053 938 1200 977">See page</th> </tr> </thead> <tbody> <tr> <td data-bbox="514 977 1053 1016">About the Analysis Parameters</td> <td data-bbox="1053 977 1200 1016">5-17</td> </tr> <tr> <td data-bbox="514 1016 1053 1058">Using Analysis Parameter Files</td> <td data-bbox="1053 1016 1200 1058">5-24</td> </tr> </tbody> </table>	Topic	See page	About the Analysis Parameters	5-17	Using Analysis Parameter Files	5-24
Topic	See page						
About the Analysis Parameters	5-17						
Using Analysis Parameter Files	5-24						
6	<p>Double-click the size standard text field or the analysis parameters text field to edit the size standard or the analysis parameters.</p> <p>For more information, see:</p> <table border="1" data-bbox="514 1218 1200 1338"> <thead> <tr> <th data-bbox="514 1218 1053 1256">Topic</th> <th data-bbox="1053 1218 1200 1256">See page</th> </tr> </thead> <tbody> <tr> <td data-bbox="514 1256 1053 1295">Editing an Existing Standard</td> <td data-bbox="1053 1256 1200 1295">5-37</td> </tr> <tr> <td data-bbox="514 1295 1053 1338">Changing an Existing Analysis Parameters File</td> <td data-bbox="1053 1295 1200 1338">5-27</td> </tr> </tbody> </table>	Topic	See page	Editing an Existing Standard	5-37	Changing an Existing Analysis Parameters File	5-27
Topic	See page						
Editing an Existing Standard	5-37						
Changing an Existing Analysis Parameters File	5-27						

No.	Description
7	A notation appears in this Information Display field when you move the cursor over a Sample file name or over a dye color field. For information on how to customize this field, see “Displaying Sample and Dye Information” on page 5-12.

Customizing the Display


The following table lists how you can customize the display by changing the settings:

Note These preferences also apply to the Results Control window. See Chapter 7, “Evaluating Analysis Results.”

To change...	Choose...	For more information
information displayed in the information display field	Project Options from the Settings menu and Sample Info Display from the submenu.	“Displaying Sample and Dye Information” on page 5-12.
the sorting of sample files	Project Options from the Settings menu and Sample File Sorting from the submenu.	“Setting Sample File Sort Order” on page 5-14.
dye Indicator code dye color	Preferences from the Setting menu and Dye Indicators from the submenu.	“Setting Dye Indicator Preferences” on page 5-15.

Using the Analysis Control Window

To use the Analysis Control window:

To...	Then...	Result
select all the samples	click the upper-left cell. Click here 	All the columns in the Analysis Control window are selected.
select all of one dye color	click the column heading for that dye color.	The column is highlighted for the color selected.

To use the Analysis Control window: *(continued)*

To...	Then...	Result
select all four dyes for one sample	click the row number.	All the colors in the row are selected.
change all the standards in the size standards column to the same setting	a. Click the arrow in the column heading. b. Choose a file from the pop-up menu.	The same size standard is displayed for all the samples.
change a size standard in one of the rows	a. Click the arrow in the row. b. Choose a file from the pop-up menu.	The size standard for the selected row changes.
change all the parameters in the parameters column to the same setting	a. Click the arrow in the row heading. b. Choose a file from the pop-up menu.	The same parameter is displayed for all the samples.
change a parameter in one of the rows	a. Click the arrow in the row. b. Choose a file from the pop-up menu.	The parameter for the selected row changes.
apply a choice to selected fields in the size standards or parameters column	a. Click the row in the column containing the information you want to apply and drag down. b. Choose Fill Down from the Edit menu.	The value in the selected rows changes to the value in the first row selected.

Analyzing Sample Files: Using the Analysis Control Window

Introduction This section describes using the Analysis Control window to perform the following tasks:

Topic	See page
Accessing Sample Files	5-6
Analyzing Sample Files	5-7
Specifying the Format for Printed Results	5-9
Displaying Size Standards and Analysis Parameters	5-11
Displaying Sample and Dye Information	5-12
Setting Sample File Sort Order	5-14
Setting Dye Indicator Preferences	5-15

Accessing Sample Files There are two ways to access Sample files contained in a project from the Analysis Control window.

Sample files that are dimmed could not be found by the project.

You can...	Then...						
double-click a Sample file name	<table border="1"> <thead> <tr> <th>If the Sample file is...</th> <th>Then that...</th> </tr> </thead> <tbody> <tr> <td>open</td> <td>Sample file window becomes active.</td> </tr> <tr> <td>not open</td> <td>Sample file opens to its Sample Results view.</td> </tr> </tbody> </table>	If the Sample file is...	Then that...	open	Sample file window becomes active.	not open	Sample file opens to its Sample Results view.
	If the Sample file is...	Then that...					
	open	Sample file window becomes active.					
not open	Sample file opens to its Sample Results view.						
select a Sample file and choose one of the five display modes from the Sample menu	the Sample file window appears in the display mode selected.						

Analyzing Sample Files

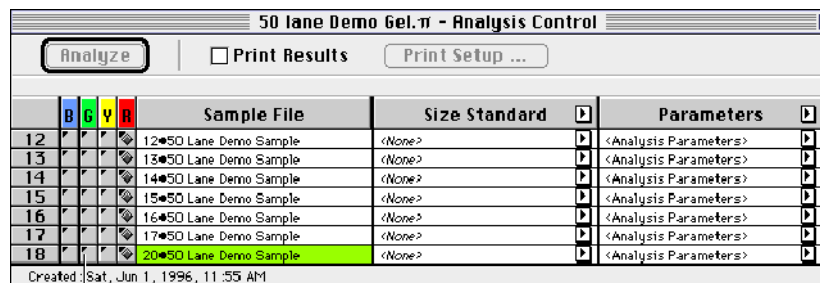
The Analysis Control window allows you to analyze multiple samples easily. You choose dyes to analyze, dye standard, size standard, and analysis parameters for each Sample file, and then analyze the Sample files using these settings. You can also install a new matrix to the Sample files.

To analyze Sample files:

Step	Action								
1	Click the dye color fields for each sample you want to analyze as follows:								
	<table border="1"> <thead> <tr> <th>To...</th> <th>Click the...</th> </tr> </thead> <tbody> <tr> <td>select a dye for all samples</td> <td>colored column header for that dye.</td> </tr> <tr> <td>select all dyes for a single Sample file</td> <td>index number at the left end of the row in which the Sample file appears.</td> </tr> <tr> <td>all dyes for all samples</td> <td>area above the row index numbers.</td> </tr> </tbody> </table>	To...	Click the...	select a dye for all samples	colored column header for that dye.	select all dyes for a single Sample file	index number at the left end of the row in which the Sample file appears.	all dyes for all samples	area above the row index numbers.
	To...	Click the...							
	select a dye for all samples	colored column header for that dye.							
select all dyes for a single Sample file	index number at the left end of the row in which the Sample file appears.								
all dyes for all samples	area above the row index numbers.								
select a dye for all samples	colored column header for that dye.								
select all dyes for a single Sample file	index number at the left end of the row in which the Sample file appears.								
all dyes for all samples	area above the row index numbers.								
2	Identify the sample containing the standard as follows:								
	<table border="1"> <thead> <tr> <th>To...</th> <th>⌘-click the...</th> </tr> </thead> <tbody> <tr> <td>identify each sample that contains a standard</td> <td>colored field that represents the standard.</td> </tr> <tr> <td>select the same dye as the standard for all samples</td> <td>colored column heading for that dye.</td> </tr> </tbody> </table>	To...	⌘-click the...	identify each sample that contains a standard	colored field that represents the standard.	select the same dye as the standard for all samples	colored column heading for that dye.		
	To...	⌘-click the...							
	identify each sample that contains a standard	colored field that represents the standard.							
select the same dye as the standard for all samples	colored column heading for that dye.								
identify each sample that contains a standard	colored field that represents the standard.								
select the same dye as the standard for all samples	colored column heading for that dye.								
A gray diamond appears in the field to identify the dye color as the standard.									
3	Choose a defined size standard setting from the pop-up menu in the Size Standard column as follows:								
	<table border="1"> <thead> <tr> <th>To...</th> <th>See...</th> </tr> </thead> <tbody> <tr> <td>define a new size standard</td> <td>"Using Size Standards" on page 5-36.</td> </tr> <tr> <td>edit a size standard</td> <td>"Editing an Existing Standard" on page 5-37.</td> </tr> <tr> <td>change the available size standards in the pop-up menu</td> <td>"Defining Folder Locations" on page 5-41.</td> </tr> </tbody> </table>	To...	See...	define a new size standard	"Using Size Standards" on page 5-36.	edit a size standard	"Editing an Existing Standard" on page 5-37.	change the available size standards in the pop-up menu	"Defining Folder Locations" on page 5-41.
	To...	See...							
	define a new size standard	"Using Size Standards" on page 5-36.							
edit a size standard	"Editing an Existing Standard" on page 5-37.								
change the available size standards in the pop-up menu	"Defining Folder Locations" on page 5-41.								
define a new size standard	"Using Size Standards" on page 5-36.								
edit a size standard	"Editing an Existing Standard" on page 5-37.								
change the available size standards in the pop-up menu	"Defining Folder Locations" on page 5-41.								

To analyze Sample files: (continued)

Step	Action						
4	To install a new matrix, select a set or all of the samples in the Sample File column and choose Install New Matrix from the Sample Menu.						
5	Choose a parameters setting from the pop-up menu in the Parameters column as follows:						
	<table border="1"> <thead> <tr> <th>To...</th> <th>See...</th> </tr> </thead> <tbody> <tr> <td>use the default analysis parameters to specify different parameters</td> <td>"Defining Analysis Parameters" on page 5-18.</td> </tr> <tr> <td>change the available parameters in the pop-up menu</td> <td>"Defining Folder Locations" on page 5-41.</td> </tr> </tbody> </table>	To...	See...	use the default analysis parameters to specify different parameters	"Defining Analysis Parameters" on page 5-18.	change the available parameters in the pop-up menu	"Defining Folder Locations" on page 5-41.
	To...	See...					
use the default analysis parameters to specify different parameters	"Defining Analysis Parameters" on page 5-18.						
change the available parameters in the pop-up menu	"Defining Folder Locations" on page 5-41.						
6	Select the Print Results checkbox to print the results automatically. For information on print set-up, see "Specifying the Format for Printed Results" on page 5-9.						
7	Click Analyze.						
	<table border="1"> <thead> <tr> <th>If a field...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>contains a small triangle (see the figure below)</td> <td>you have analyzed the samples using the currently selected size standard and analysis parameters.</td> </tr> <tr> <td>does not contain a triangle</td> <td>the sample has not yet been analyzed with the current settings.</td> </tr> </tbody> </table>	If a field...	Then...	contains a small triangle (see the figure below)	you have analyzed the samples using the currently selected size standard and analysis parameters.	does not contain a triangle	the sample has not yet been analyzed with the current settings.
	If a field...	Then...					
contains a small triangle (see the figure below)	you have analyzed the samples using the currently selected size standard and analysis parameters.						
does not contain a triangle	the sample has not yet been analyzed with the current settings.						



Triangles indicate samples were analyzed with the current setting

To analyze Sample files: *(continued)*

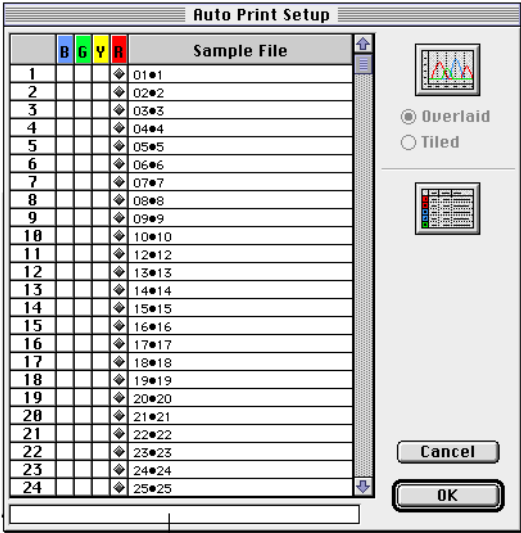
Step	Action
8	<p>The triangle disappears if you</p> <ul style="list-style-type: none">◆ Select a different size standard or analysis parameters file for a dye/sample.◆ Modify the currently selected size standard or analysis parameters file (or the analysis parameters stored as program preferences).◆ Specify a different dye color as the standard (move the diamond indicator to another dye/sample).◆ Assign a new matrix to the Sample file.
9	<p>To verify the results, refer to “Process of Verifying Results” on page 7-46.</p>

Specifying the Format for Printed Results

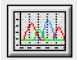

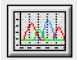

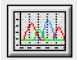

To specify the format for printed results:

Step	Action
1	<p>In the Analysis Control window, select the checkbox labeled Print Results.</p> <p>When this checkbox is selected, the Print Setup button becomes active.</p>
2	<p>Click the Print Setup button.</p> <p>The Auto Print Setup dialog box appears (see below).</p> <p>All the dyes selected for analysis are also selected for printing.</p>

To specify the format for printed results: *(continued)*

Step	Action
	<p data-bbox="481 185 733 211">Sample Dye color fields</p>  <p data-bbox="481 782 830 808">Sample Information Display field</p>
3	Moving the cursor over a Sample file name or over a dye color field, a notation appears in the Sample Information Display field.
4	Click the sample dye color fields to specify any sample you do not want to print (see the above figure).

To specify the format for printed results: *(continued)*

Step	Action						
5	Choose the format by clicking either or both of the buttons at the right of the window.						
	<table border="1"> <thead> <tr> <th>Click this button...</th> <th>To print...</th> </tr> </thead> <tbody> <tr> <td>  </td> <td> electropherograms for the samples and dyes selected for analysis. Select the appropriate radio button to specify whether the electropherograms for the four dyes appear: <ul style="list-style-type: none"> ◆ Together in one panel (overlaid), or in ◆ Separate panels (tiled). </td> </tr> <tr> <td>  </td> <td>tabular data.</td> </tr> </tbody> </table>	Click this button...	To print...		electropherograms for the samples and dyes selected for analysis. Select the appropriate radio button to specify whether the electropherograms for the four dyes appear: <ul style="list-style-type: none"> ◆ Together in one panel (overlaid), or in ◆ Separate panels (tiled). 		tabular data.
	Click this button...	To print...					
	electropherograms for the samples and dyes selected for analysis. Select the appropriate radio button to specify whether the electropherograms for the four dyes appear: <ul style="list-style-type: none"> ◆ Together in one panel (overlaid), or in ◆ Separate panels (tiled). 						
	tabular data.						
6	Click OK.						

Displaying Size Standards and Analysis Parameters

Use the Analysis Control window to open, review, or change Size Standard and analysis parameters.

To display these files:

You can...	Then...	For more information see...
double click the field containing the size standard or the analysis parameters file	the Size Standard or Analysis Parameter window appears.	<ul style="list-style-type: none"> ◆ "Using Analysis Parameter Files" on page 5-24. ◆ "Defining the Size Standard" on page 5-31.

Displaying Sample and Dye Information

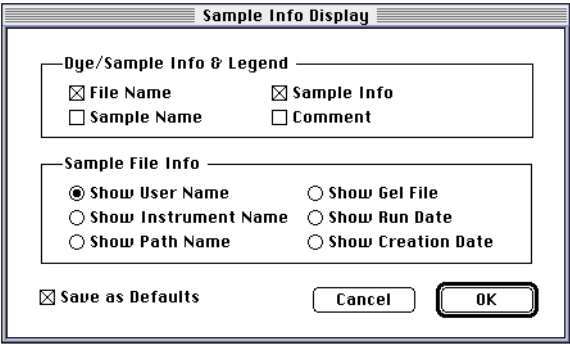
How to Display Sample and Dye Information

Move the cursor over a dye color field or over a Sample file name field to display information about the samples and dyes.

How to Specify the Information Displayed

The following procedure describes how to specify the information displayed when moving the cursor over a dye color field or a Sample file name field.

To specify the information displayed:

Step	Action										
1	<p>Choose Project Options from the Settings menu and Sample Info Display from the submenu.</p> <p>The Sample Info Display dialog box appears.</p> 										
2	<p>Select the checkboxes in the Dye/Sample Info & Legend section to control what appears when you move the cursor over the dye color fields in the Control windows.</p> <p>The following table describes the checkboxes:</p> <table border="1" data-bbox="467 1188 1189 1445"> <thead> <tr> <th>If you select this checkbox...</th> <th>This appears...</th> </tr> </thead> <tbody> <tr> <td>File Name</td> <td>Sample file name.</td> </tr> <tr> <td>Sample Name</td> <td>Name of the sample file from the Sample file.</td> </tr> <tr> <td>Sample Info</td> <td>Sample information from the Sample file.</td> </tr> <tr> <td>Comment</td> <td>Comment from the Sample file.</td> </tr> </tbody> </table>	If you select this checkbox...	This appears...	File Name	Sample file name.	Sample Name	Name of the sample file from the Sample file.	Sample Info	Sample information from the Sample file.	Comment	Comment from the Sample file.
If you select this checkbox...	This appears...										
File Name	Sample file name.										
Sample Name	Name of the sample file from the Sample file.										
Sample Info	Sample information from the Sample file.										
Comment	Comment from the Sample file.										

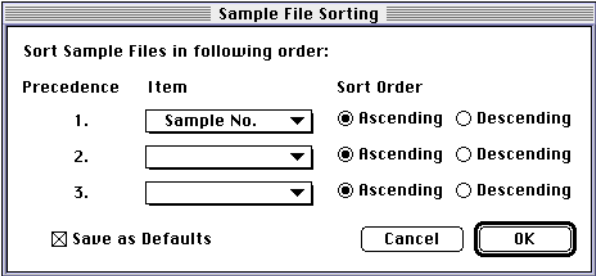
To specify the information displayed: *(continued)*

Step	Action														
3	<p>In the Sample File Info heading, select the checkboxes for the information that you want to display when you move the cursor over the Sample File name field.</p> <p>The following table describes the radio buttons:</p> <table border="1" data-bbox="510 349 1229 803"> <thead> <tr> <th data-bbox="517 357 873 389">Select this radio button...</th> <th data-bbox="873 357 1223 389">To display...</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 389 873 462">Show User Name</td> <td data-bbox="873 389 1223 462">user name from the Sample file.</td> </tr> <tr> <td data-bbox="517 462 873 527">Show Instrument Name</td> <td data-bbox="873 462 1223 527">instrument name from the Sample file.</td> </tr> <tr> <td data-bbox="517 527 873 592">Show Folder Name</td> <td data-bbox="873 527 1223 592">path and name of the folder where the file is located.</td> </tr> <tr> <td data-bbox="517 592 873 657">Show Gel File</td> <td data-bbox="873 592 1223 657">name of gel file from the Sample file.</td> </tr> <tr> <td data-bbox="517 657 873 722">Show Run Date</td> <td data-bbox="873 657 1223 722">run date and start time from the Sample file.</td> </tr> <tr> <td data-bbox="517 722 873 795">Show Creation Date</td> <td data-bbox="873 722 1223 795">date and time the Sample file was created.</td> </tr> </tbody> </table>	Select this radio button...	To display...	Show User Name	user name from the Sample file.	Show Instrument Name	instrument name from the Sample file.	Show Folder Name	path and name of the folder where the file is located.	Show Gel File	name of gel file from the Sample file.	Show Run Date	run date and start time from the Sample file.	Show Creation Date	date and time the Sample file was created.
Select this radio button...	To display...														
Show User Name	user name from the Sample file.														
Show Instrument Name	instrument name from the Sample file.														
Show Folder Name	path and name of the folder where the file is located.														
Show Gel File	name of gel file from the Sample file.														
Show Run Date	run date and start time from the Sample file.														
Show Creation Date	date and time the Sample file was created.														
4	Select the checkbox labeled Save as Defaults to have the options you choose saved as the default settings.														
5	Click OK.														

Setting Sample File Sort Order

Use this option to specify how Sample files are sorted using three criteria. If no sorting option is specified, then the program sorts by Sample number.

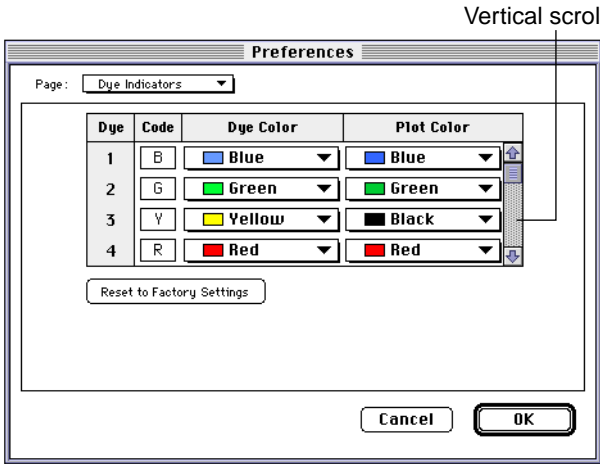
To set Sample file sort order:

Step	Action
1	<p>Choose Project Options from the Settings menu and Sample File Sorting from the submenu. The Sample File Sorting dialog box appears.</p> 
2	<p>Choose from the following items from the pop-up menus:</p> <ul style="list-style-type: none"> ◆ File Name. ◆ Directory. ◆ Gel Name. ◆ Sample Number. ◆ User Name. ◆ Instrument Name. ◆ Run Data. ◆ Creation Date. ◆ As Added (sorts the files in the order that they were added). <p>The precedence indicates the sorting level.</p>
3	Select a radio button for each sort to indicate whether ascending or descending order.
4	Select the checkbox labeled Save as Defaults to have the options you choose saved as the default settings.
5	Click OK.

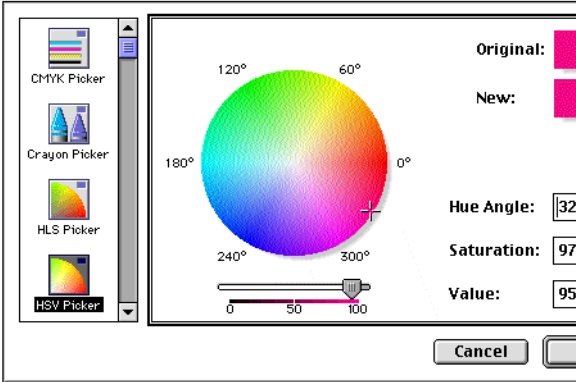
Setting Dye Indicator Preferences

The following procedure describes how to change the defaults that determine what dye colors appear on the screen and on printed results. Setting default dye and plot colors sets the colors used for both the Control windows and the Results displays.

To set default dye and plot colors:

Step	Action						
1	<p>Choose Preferences from the Settings menu and Dye Indicators from the submenu. The Preferences window appears.</p> <p>Note If the Preferences window is already displayed, choose Dye Indicators from the pop-up menu.</p> 						
2	<p>The following table describes the Dye Color and Plot Color columns:</p> <table border="1"> <thead> <tr> <th>Item</th> <th>Description</th> </tr> </thead> <tbody> <tr> <td>Dye Color column</td> <td>Shows the colors that represent the dyes in the Control window lists and the gel display of ABI™ 373 and ABI PRISM® 377 data. The dye color is also identified in the left color legend in the Results display.</td> </tr> <tr> <td>Plot Color column</td> <td>Shows the colors used for plotting the data in the electropherograms.</td> </tr> </tbody> </table>	Item	Description	Dye Color column	Shows the colors that represent the dyes in the Control window lists and the gel display of ABI™ 373 and ABI PRISM® 377 data. The dye color is also identified in the left color legend in the Results display.	Plot Color column	Shows the colors used for plotting the data in the electropherograms.
Item	Description						
Dye Color column	Shows the colors that represent the dyes in the Control window lists and the gel display of ABI™ 373 and ABI PRISM® 377 data. The dye color is also identified in the left color legend in the Results display.						
Plot Color column	Shows the colors used for plotting the data in the electropherograms.						
3	Use the vertical scroll bar to change the dye color and plot color for a fifth dye.						

To set default dye and plot colors: *(continued)*

Step	Action						
4	To change a code, type a different character in the appropriate entry field in the Code column.						
5	<p>To change a color, select a new color from the pop-up menu.</p> <p>If you choose Other from the pop-up menu, a color picker appears.</p>  <p>Using the color picker:</p> <table border="1" data-bbox="467 841 1192 1203"> <thead> <tr> <th data-bbox="467 841 655 878">To use the...</th> <th data-bbox="655 841 1192 878">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="467 878 655 1052">scrolling fields</td> <td data-bbox="655 878 1192 1052"> <p>click one of the up or down arrows and hold down the mouse button.</p> <p>The new color appears in the New box and the Original box retains the original color, until clicking OK to accept the new color.</p> </td> </tr> <tr> <td data-bbox="467 1052 655 1203">color wheel</td> <td data-bbox="655 1052 1192 1203"> <ol style="list-style-type: none"> a. Click the desired color in the wheel. The new color appears in the New box. b. Move the insertion point and click again until the desired color appears. </td> </tr> </tbody> </table>	To use the...	Then...	scrolling fields	<p>click one of the up or down arrows and hold down the mouse button.</p> <p>The new color appears in the New box and the Original box retains the original color, until clicking OK to accept the new color.</p>	color wheel	<ol style="list-style-type: none"> a. Click the desired color in the wheel. The new color appears in the New box. b. Move the insertion point and click again until the desired color appears.
To use the...	Then...						
scrolling fields	<p>click one of the up or down arrows and hold down the mouse button.</p> <p>The new color appears in the New box and the Original box retains the original color, until clicking OK to accept the new color.</p>						
color wheel	<ol style="list-style-type: none"> a. Click the desired color in the wheel. The new color appears in the New box. b. Move the insertion point and click again until the desired color appears. 						
6	Click OK when finished changing dye indicator preferences.						

About the Analysis Parameters

Introduction The GeneScan Analysis Software has default analysis parameters that are stored as preferences. You can set and use these default analysis parameters, or create analysis parameters files to use with specific protocols.

Parameters Used During Automatic Analysis During automatic analysis, the GeneScan Analysis Software uses the information from the data collection Injection List or Run Sheet window to determine which parameters to use.

If you do not specify the information in the Data Collection software, the GeneScan Analysis Software uses the parameters set in the Auto-Analysis Defaults (see step 2 on page 4-6).

Specifying Where Files are Stored Choose Preferences from the Settings menu and Folder Locations from the submenu to specify the folders in which the software automatically stores the analysis parameters and size standards files.

For more information, see “Defining Folder Locations” on page 5-41.

Defining Analysis Parameters

Choose Analysis Parameters from the Settings menu to define the Analysis Parameters.

The Analysis Parameters dialog box appears.

Analysis Parameters

Analysis Range

Full Range
 This Range (Data Points)
Start:
Stop:

Size Call Range

All Sizes
 This Range (Base Pairs)
Min:
Max:

Data Processing

Baseline
 MultiComponent

Smooth Options

None
 Light
 Heavy

Peak Detection

Peak Amplitude Thresholds

B: Y:
G: R:

Min. Peak Half Width: Pts

Size Calling Method

2nd Order Least Squares
 3rd Order Least Squares
 Cubic Spline Interpolation
 Local Southern Method
 Global Southern Method

Split Peak Correction

None
 GENESCAN 2500
 LeftMost Peak
 RightMost Peak

Correction Limit: Data Pts

Cancel OK

Six Analysis Parameters

There are six Analysis parameters:

Topic	See page
Analysis Range Parameter Options	5-19
Data Processing Parameter Options	5-19
Peak Detection Parameter Options	5-20
Size Call Range Parameter Options	5-21
Size Calling Method Parameter Options	5-21
Split Peak Correction Parameter Options	5-22

Analysis Range Parameter Options

The following are the Analysis Range parameter options:

Item	Description
Full Range radio button	Use to analyze all the data collected on the genetic analysis instrument for each sample.
This Range (Data Points) radio button	<p>Enter Start and Stop data point numbers in the entry fields in order to specify only a limited range analyzed for each sample.</p> <p>This affects what is displayed in the results display. Normally, set the analysis range to start after the primer peak.</p> <p>Note Sample files generated from ABI™ 373 or ABI PRISM® 377 may have already removed the primer peak. This was done by setting the scan range for gel image generation to exclude the primer peak.</p>

Data Processing Parameter Options

About the Data Processing Parameter Options

The Data Processing parameter options specify how the raw data is processed before peak detection and size calling.

Data Processing Parameter Options Described

Item	Description
Baseline checkbox	<p>Use to automatically adjust the baselines of all detected dye colors to the same level for a better comparison of relative signal intensity.</p> <p>When the GeneScan Analysis Software data is sent to the Macintosh® computer, during data collection, the baseline for each dye is at a different level.</p>
Multicomponent checkbox	<p>Use to specify that the GeneScan Analysis Software apply a predefined matrix to adjust for spectral overlap when it performs analysis.</p> <p>Although the dyes used to label DNA fluoresce at different wavelengths, the spectra overlap to some extent. Create a matrix file containing a mathematical matrix to correct for this overlap.</p> <p>For a description of matrix files and how to create them, see Chapter 6, "Making a Matrix File."</p>

Item	Description								
Smooth Options	Use to help reduce the number of false peaks detected by the GeneScan Analysis Software. You have the following options:								
	<table border="1"> <thead> <tr> <th>Select...</th> <th>To...</th> </tr> </thead> <tbody> <tr> <td>None</td> <td>apply no smoothing. Select this option if the data has very sharp, narrow peaks of interest.</td> </tr> <tr> <td>Light</td> <td>provide the best results for normal data.</td> </tr> <tr> <td>Heavy</td> <td>apply to data from slower runs that has very broad peaks. Choosing this option might reduce peak size or eliminate narrow peaks.</td> </tr> </tbody> </table>	Select...	To...	None	apply no smoothing. Select this option if the data has very sharp, narrow peaks of interest.	Light	provide the best results for normal data.	Heavy	apply to data from slower runs that has very broad peaks. Choosing this option might reduce peak size or eliminate narrow peaks.
	Select...	To...							
	None	apply no smoothing. Select this option if the data has very sharp, narrow peaks of interest.							
Light	provide the best results for normal data.								
Heavy	apply to data from slower runs that has very broad peaks. Choosing this option might reduce peak size or eliminate narrow peaks.								

Peak Detection Parameter Options

About the Peak Detection Parameter Options

Use the Peak Detection parameter options to specify the minimum peak height to be detected for analysis. This, in turn, controls the number of peaks analyzed. Peaks falling below the parameters specified appear in the electropherogram, but are not analyzed, and no values appear for them in the tabular data.

Peak Detection Parameter Options Described

Item	Description	For example
Dye Amplitude Threshold	Set the dye amplitude threshold at a level that allows the software to detect peaks, but eliminate noise. For each dye, the GeneScan Analysis Software detects peaks above the threshold entered in the entry field.	If you leave the default value of 50, peaks with amplitude above 50 are analyzed and appear in the tabular data. Lower amplitude peaks still appear in the electropherogram, but are not analyzed and do not appear in the tabular data.

Item	Description	For example
Minimum Peak Half Width	<p>Defines what constitutes a peak.</p> <p>Use to specify the smallest half peak width for peak detection.</p> <p>The range is from 2 - 99.</p> <p>A typical number might be 3 for microsattelites, or 10 for SSCPs.</p>	<p>If this number is large, the software ignores noise spikes.</p> <p>If the peaks in the data are narrow, set the value to a low number.</p> <p>Experiment with this value to determine the best number for the data.</p>

Size Call Range Parameter Options

About the Size Call Range Parameter Options

Use the Size Call Range parameter options to specify the range of size fragment (in base pairs) to be included in the peak tabular data.

Size Call Range Parameter Options Described

Item	Description
All Sizes radio button	All detected fragments appear in the tabular data.
This Range (Base Pairs) radio button	Define lower (minimum) and upper (maximum) limits of the peaks to include in the tabular data.

Size Calling Method Parameter Options

About Size Calling Method Parameter Options

Click a radio button to select the desired size calling method. The GeneScan Analysis Software uses these methods to determine the molecular length of an unknown fragment.

Size Calling Method Parameter Options Described

Item	Description
2nd Order Least Squares and 3rd Order Least Squares	<p>Both Least Squares methods use regression analysis to build a best-fit size calling curve.</p> <p>For more information, see "Least Square Method" on page C-2.</p>

Item	Description
Cubic Spline Interpolation	Forces the sizing curve through all the known points of the selected GeneScan size standard. For more information, see “Cubic Spline Interpolation Method” on page C-4.
Local Southern method	Determines the sizes of fragments by using the reciprocal relationship between fragment length and mobility. For more information, see “Local Southern Method” on page C-5.
Global Southern method	Similar to the Least Squares method in that it compensates for standard fragments that may run anomalously. For more information, see “Global Southern Method” on page C-7.

Split Peak Correction Parameter Options

About the Split Peak Feature

Under denaturing conditions, certain fragments in the GeneScan 2500 standard appear as doublets, or split peaks. This standard has labels on both stands of the DNA. Under poor denaturing conditions you see split peaks. One of the two fragments typically has normal mobility, while the other does not. The Split Peak Correction feature allows the software to correctly call each of the splits.

Split Peak Parameter Options Described

If you select split peak correction, you will also need to verify or change the correction limits. After you decide on a split peak correction method, use the same method for all projects to keep size calling consistent.

Split Peak Correction parameter options:

Item	Description
None	No correction for doublets.
GeneScan 2500	Makes the following peak size assignments for GeneScan 2500: <ul style="list-style-type: none"> ◆ The right peak for all fragments 222, 233, 238, 286, and 490. ◆ The left peak for all other splits.
LeftMost Peak	Chooses the left peak for every doublet.

Split Peak Correction parameter options: *(continued)*

Item	Description
RightMost Peak	Chooses the right peak for every doublet.
Correction Limit	Set a correction limit if correcting for doublets. Set this value slightly larger than the largest split observed. This value (set in scan lines or data points) specifies the maximum width of split that should be corrected (the difference in scan numbers or data points of the positions of the two peaks).

Using Analysis Parameter Files

In This Section This section contains the following topics:

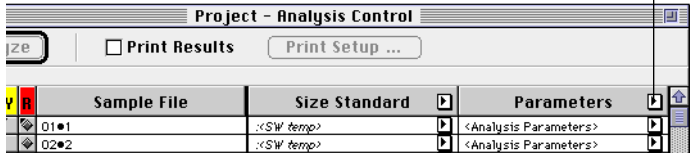
Topic	See page
Analyzing Samples Using the Same Analysis Parameters	5-24
Selecting Different Parameters for Analysis Samples	5-25
Displaying Default Parameters	5-26
Creating Custom Analysis Parameter Files	5-26
Changing an Existing Analysis Parameters File	5-27
Deleting Custom Analysis Parameters	5-28

Analyzing Samples Using the Same Analysis Parameters

To analyze samples using the same analysis parameters:

Step	Action
1	If the Analysis Control window is not displayed, then choose Analysis Control (⌘ 1) from the Windows menu.
2	Click the arrow in the Parameters column heading and choose parameters from the pop-up menu. Your menu choice applies to all fields in the column.

Pop-up menu



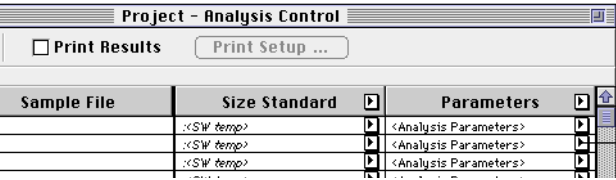
The screenshot shows the 'Project - Analysis Control' window. It contains a table with the following data:

Sample File	Size Standard	Parameters
01#1	:S\W temp	<Analysis Parameters>
02#2	:S\W temp	<Analysis Parameters>

Step	Action								
3	<p>The pop-up menu contains the following options:</p> <table border="1"> <thead> <tr> <th>Item</th> <th>Description</th> </tr> </thead> <tbody> <tr> <td><Analysis Parameters></td> <td>Applies the parameters that are stored as preferences in the software.</td> </tr> <tr> <td><Collection Settings></td> <td>Applies the analysis parameters file specified in the data collection program, which is embedded in the Sample file.</td> </tr> <tr> <td>Custom parameters that are listed at the bottom of the menu</td> <td> <p>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p>For information on setting the Folder Locations preferences, see “Specifying File Locations” on page 5-42.</p> </td> </tr> </tbody> </table>	Item	Description	<Analysis Parameters>	Applies the parameters that are stored as preferences in the software.	<Collection Settings>	Applies the analysis parameters file specified in the data collection program, which is embedded in the Sample file.	Custom parameters that are listed at the bottom of the menu	<p>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p>For information on setting the Folder Locations preferences, see “Specifying File Locations” on page 5-42.</p>
Item	Description								
<Analysis Parameters>	Applies the parameters that are stored as preferences in the software.								
<Collection Settings>	Applies the analysis parameters file specified in the data collection program, which is embedded in the Sample file.								
Custom parameters that are listed at the bottom of the menu	<p>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p>For information on setting the Folder Locations preferences, see “Specifying File Locations” on page 5-42.</p>								

Selecting Different Parameters for Analysis Samples

To apply separate analysis parameters to selected samples:

Step	Action
1	If the Analysis Control window is not displayed, then choose Analysis Control (⌘ 1) from the Windows menu.
2	<p>Click the arrow in the Parameters column for the sample that you want to change the parameter settings.</p> <p>A pop-up menu appears.</p>  <p>The screenshot shows the 'Project - Analysis Control' window. At the top, there are buttons for 'Print Results' and 'Print Setup ...'. Below is a table with three columns: 'Sample File', 'Size Standard', and 'Parameters'. The 'Parameters' column has a dropdown arrow for each row. A 'Pop-up menu' is shown over the first row's dropdown arrow, displaying options: '<Analysis Parameters>', '<Analysis Parameters>', and '<Analysis Parameters>'. The 'Sample File' column contains entries like ':<SW temp>'. The 'Size Standard' column contains a dropdown arrow. The 'Parameters' column contains a dropdown arrow and a small icon.</p>

Step	Action						
3	The pop-up menu contains the following options:						
	<table border="1"> <thead> <tr> <th>Choose...</th> <th>To...</th> </tr> </thead> <tbody> <tr> <td><Analysis Parameters></td> <td>apply the parameters that are stored as preferences in the software.</td> </tr> <tr> <td>Define New</td> <td>display the Analysis Control dialog box. For information on completing the fields, see "About the Analysis Parameters" on page 5-17.</td> </tr> </tbody> </table>	Choose...	To...	<Analysis Parameters>	apply the parameters that are stored as preferences in the software.	Define New	display the Analysis Control dialog box. For information on completing the fields, see "About the Analysis Parameters" on page 5-17.
	Choose...	To...					
<Analysis Parameters>	apply the parameters that are stored as preferences in the software.						
Define New	display the Analysis Control dialog box. For information on completing the fields, see "About the Analysis Parameters" on page 5-17.						
4	Repeat step 2 and step 3 for each sample. Note You can also use the Macintosh computer Cut, Copy, and Paste commands from the Edit menu.						

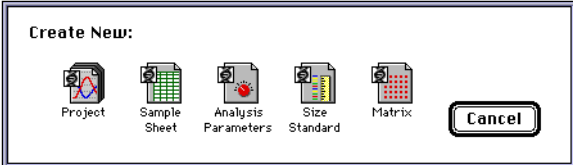
Displaying Default Parameters

There are two ways to display the default parameters:

You can...	Then...
choose Analysis Parameters from the Settings menu.	Analysis Parameters dialog box appears with the default parameters.
double-click <Analysis Parameters> from within the Analysis Control window.	For information on defining the Analysis Parameters, see page 5-18.

Creating Custom Analysis Parameter Files

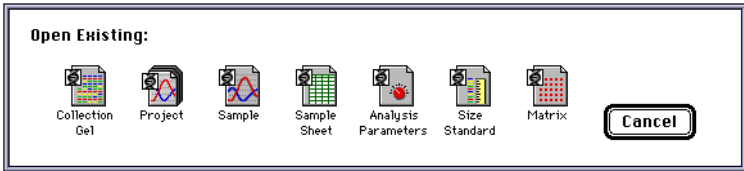
To create a custom analysis parameter file:

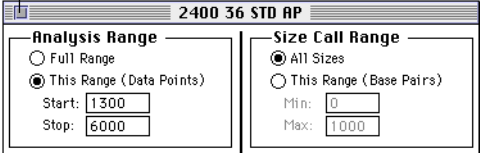
Step	Action
1	<p>Choose New from the File menu.</p> <p>The Create New dialog box appears.</p> 

Step	Action
2	Click the Analysis Parameters icon. The Analysis Parameters dialog box appears.
3	Change the parameters as necessary. For information on the specific options, see “About the Analysis Parameters” on page 5-17.
4	Choose Save (⌘ S) from the File menu. A dialog box appears.
5	Enter a descriptive name and click Save. Note Unless you choose a different folder location, the GeneScan Analysis Software stores the file in the folder specified as the folder location for Analysis Parameters. For more information, see “Defining Folder Locations” on page 5-41. The file now appears in the pop-up menu for analysis parameters in the GeneScan Analysis Control window. You can also select the file in the Data Collection software for automatic analysis.

Changing an Existing Analysis Parameters File

To change an existing analysis parameter file:

Step	Action
1	Choose Open (⌘ O) from the File menu. Note You can also double-click the file name in the Parameters column of the Analysis Control window. The Open Existing dialog box appears.
	
2	Click the Analysis Parameters icon. A directory dialog box appears.
3	Select a file that you want to change and click Open.

Step	Action
4	<p>Make the changes and close the window by clicking the close box.</p> <p>Close box</p> 

Deleting Custom Analysis Parameters

To delete a custom analysis parameters file:

Step	Action
1	<p>Find the file in the folder you specified as the location for the analysis parameters files.</p> <p>Note The custom analysis parameters file is in the folder that is normally called the GS Parameters Folder. This folder is inside the ABI PRISM GeneScan folder.</p>
2	<p>Drag the custom analysis parameters file to the trash and empty the trash.</p>

About Size Standards

What are Size Standards Size standards are specific DNA fragments of known sizes. After defining the peaks of a size standard, the GeneScan Analysis Software matches this definition to the internal size standard included with the run. The software assigns the defined size values to the appropriate peaks of the internal size standard, and uses this information with the selected size calling method to size all unknown fragments.

Advantages of Using Size Standard Running an internal size standard results in accurate and precise molecular length determination. This is because the internal size standard and the unknown fragments undergo exactly the same electrophoretic forces. The GeneScan Analysis Software can then compensate for band-shift artifacts caused by variations in the gel and in the sample from lane to lane or injection to injection.

Size Standards Provided PE Applied Biosystems provides several fluorescently labeled standards, which are described in Appendix B. You can also label and use other fragments if they better suit the fragment sizes with which you are working.

When to Define Size Standards Normally, a size standard is defined using the GeneScan Analysis Software after running the standard with samples on the instrument. The software detects peaks for a selected dye color in a selected Sample file and allows you to define the peak sizes. You can save the defined standard in a file and use it to automatically analyze other samples run with the same standard and under the same conditions.

If Split Peaks Appear Split peaks might appear in size standards in which both strands are labeled, for example GeneScan 1000 (see page B-8) and GeneScan 2500 (see page B-10). For some peaks, the two strands migrate at different rates when running under denaturing conditions, and they appear as two peaks approximately half the height of normal non-split peaks. One peak of the two runs is true to size. You should assign a size to that peak for the standard definition, and assign zero to the other peak.

For more information, see “Split Peak Correction Parameter Options” on page 5-22.

**Specifying Where
Files are Stored**

Choose Preferences from the Settings menu and Folder Locations from the submenu to specify the folders in which the software automatically stores the analysis parameters and size standards files.

For more information, see “Defining Folder Locations” on page 5-41.

Defining the Size Standard

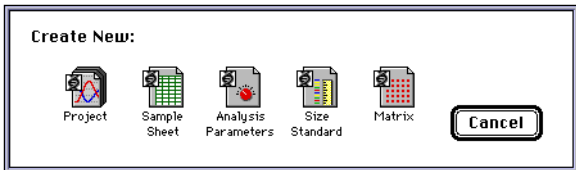
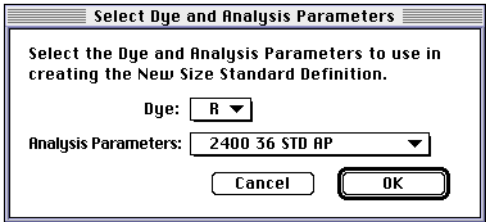
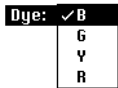
Two Ways to Define the Size Standard

There are two ways to define a new size standard:

Topic	See page
Using the New Command	5-31
Using the Analysis Control Window	5-34

Using the New Command

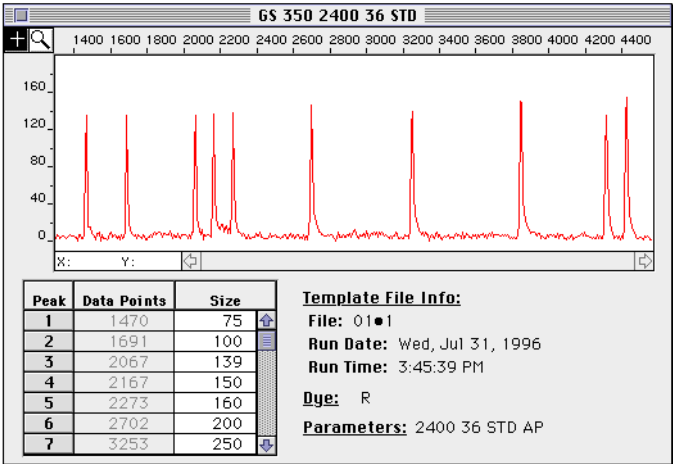
To use the New command to define a new size standard:

Step	Action
1	<p>Choose New from the File menu.</p> <p>The Create New dialog box appears.</p> 
2	<p>Click the Size Standard icon. A directory dialog box appears.</p>
3	<p>Select the Sample file that contains the dye standard you want to use as the template and click Open.</p> <p>The Select Dye and Analysis Parameters dialog box appears.</p> 
4	<p>Choose from the Dye pop-up menu the code that represents the dye label of the size standard in the selected Sample file.</p> 

To use the New command to define a new size standard: *(continued)*

Step	Action			
<p data-bbox="400 183 420 207">5</p>	<p data-bbox="471 183 1157 240">Choose from the Analysis Parameters pop-up menu the analysis parameters to use.</p> <p data-bbox="471 256 989 280">The pop-up menu contains the following options:</p>			
	<table border="1"> <thead> <tr> <th data-bbox="463 334 827 365">Item</th> <th data-bbox="827 334 1178 365">Description</th> </tr> </thead> </table>	Item	Description	
	Item	Description		
<table border="1"> <tbody> <tr> <td data-bbox="463 371 827 464"><Analysis Parameters></td> <td data-bbox="827 371 1178 464">Applies the parameters that are stored as preferences in the software.</td> </tr> <tr> <td data-bbox="463 464 827 750">custom parameters that are listed at the bottom of the menu</td> <td data-bbox="827 464 1178 750"> <p data-bbox="841 472 1173 610">These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p data-bbox="841 634 1173 740">For information on setting the Folder Locations preferences, see “Defining Folder Locations” on page 5-41.</p> </td> </tr> </tbody> </table>	<Analysis Parameters>	Applies the parameters that are stored as preferences in the software.	custom parameters that are listed at the bottom of the menu	<p data-bbox="841 472 1173 610">These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p data-bbox="841 634 1173 740">For information on setting the Folder Locations preferences, see “Defining Folder Locations” on page 5-41.</p>
<Analysis Parameters>	Applies the parameters that are stored as preferences in the software.			
custom parameters that are listed at the bottom of the menu	<p data-bbox="841 472 1173 610">These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p data-bbox="841 634 1173 740">For information on setting the Folder Locations preferences, see “Defining Folder Locations” on page 5-41.</p>			
<p data-bbox="400 773 420 797">6</p>	<p data-bbox="471 773 572 797">Click OK.</p> <p data-bbox="471 821 1177 878">A window appears (see below) showing the electropherogram and a table of peaks for the dye color and sample selected.</p> <p data-bbox="471 894 1177 951">You should be able to recognize the peak pattern of the standard in the electropherogram.</p> <p data-bbox="471 967 1123 1057">Note You can only change the peak size value in the right column of the table. You cannot change or rearrange the peak numbers.</p> <p data-bbox="471 1073 1163 1162">Note If too many peaks appear in the electropherogram or the baseline is too high, you might need to adjust the analysis parameters. See “Using Analysis Parameter Files” on page 5-24.</p> <p data-bbox="471 1179 1083 1235">The software assigns a number to each peak found in the electropherogram in order, from left to right.</p>			

To use the New command to define a new size standard: (continued)

Step	Action																								
	 <table border="1" data-bbox="548 467 801 646"> <thead> <tr> <th>Peak</th> <th>Data Points</th> <th>Size</th> </tr> </thead> <tbody> <tr> <td>1</td> <td>1470</td> <td>75</td> </tr> <tr> <td>2</td> <td>1691</td> <td>100</td> </tr> <tr> <td>3</td> <td>2067</td> <td>139</td> </tr> <tr> <td>4</td> <td>2167</td> <td>150</td> </tr> <tr> <td>5</td> <td>2273</td> <td>160</td> </tr> <tr> <td>6</td> <td>2702</td> <td>200</td> </tr> <tr> <td>7</td> <td>3253</td> <td>250</td> </tr> </tbody> </table> <p>Template File Info: File: 0101 Run Date: Wed, Jul 31, 1996 Run Time: 3:45:39 PM Dye: R Parameters: 2400 36 STD AP</p>	Peak	Data Points	Size	1	1470	75	2	1691	100	3	2067	139	4	2167	150	5	2273	160	6	2702	200	7	3253	250
Peak	Data Points	Size																							
1	1470	75																							
2	1691	100																							
3	2067	139																							
4	2167	150																							
5	2273	160																							
6	2702	200																							
7	3253	250																							
7	<p>Specify the peaks of the standards and their sizes as follows:</p> <table border="1" data-bbox="517 708 1233 1442"> <thead> <tr> <th>Step</th> <th>Action</th> </tr> </thead> <tbody> <tr> <td>a.</td> <td> <p>Click the peak you want to define either in the electropherogram or in the table.</p> <p>Use the Zoom In (⌘ Plus sign) and Zoom Out (⌘ Minus sign) commands from the View menu to zoom the electropherogram for easier viewing.</p> <table border="1" data-bbox="624 933 1228 1141"> <thead> <tr> <th>If you click a peak...</th> <th>Then the...</th> </tr> </thead> <tbody> <tr> <td>in the electropherogram</td> <td>corresponding row in the table is highlighted.</td> </tr> <tr> <td>number in the table</td> <td>corresponding peak in the electropherogram is highlighted.</td> </tr> </tbody> </table> </td> </tr> <tr> <td>b.</td> <td> <p>Type the value for the selected peak in the corresponding Size field in the table.</p> <p>Refer to Appendix B, "GeneScan Size Standards," for values and peaks patterns.</p> <p>Note Leave a zero in the Size field to ignore a peak for the standard definition.</p> </td> </tr> <tr> <td>c.</td> <td> <p>Press Return to automatically move to the next size standard peak.</p> </td> </tr> </tbody> </table>	Step	Action	a.	<p>Click the peak you want to define either in the electropherogram or in the table.</p> <p>Use the Zoom In (⌘ Plus sign) and Zoom Out (⌘ Minus sign) commands from the View menu to zoom the electropherogram for easier viewing.</p> <table border="1" data-bbox="624 933 1228 1141"> <thead> <tr> <th>If you click a peak...</th> <th>Then the...</th> </tr> </thead> <tbody> <tr> <td>in the electropherogram</td> <td>corresponding row in the table is highlighted.</td> </tr> <tr> <td>number in the table</td> <td>corresponding peak in the electropherogram is highlighted.</td> </tr> </tbody> </table>	If you click a peak...	Then the...	in the electropherogram	corresponding row in the table is highlighted.	number in the table	corresponding peak in the electropherogram is highlighted.	b.	<p>Type the value for the selected peak in the corresponding Size field in the table.</p> <p>Refer to Appendix B, "GeneScan Size Standards," for values and peaks patterns.</p> <p>Note Leave a zero in the Size field to ignore a peak for the standard definition.</p>	c.	<p>Press Return to automatically move to the next size standard peak.</p>										
Step	Action																								
a.	<p>Click the peak you want to define either in the electropherogram or in the table.</p> <p>Use the Zoom In (⌘ Plus sign) and Zoom Out (⌘ Minus sign) commands from the View menu to zoom the electropherogram for easier viewing.</p> <table border="1" data-bbox="624 933 1228 1141"> <thead> <tr> <th>If you click a peak...</th> <th>Then the...</th> </tr> </thead> <tbody> <tr> <td>in the electropherogram</td> <td>corresponding row in the table is highlighted.</td> </tr> <tr> <td>number in the table</td> <td>corresponding peak in the electropherogram is highlighted.</td> </tr> </tbody> </table>	If you click a peak...	Then the...	in the electropherogram	corresponding row in the table is highlighted.	number in the table	corresponding peak in the electropherogram is highlighted.																		
If you click a peak...	Then the...																								
in the electropherogram	corresponding row in the table is highlighted.																								
number in the table	corresponding peak in the electropherogram is highlighted.																								
b.	<p>Type the value for the selected peak in the corresponding Size field in the table.</p> <p>Refer to Appendix B, "GeneScan Size Standards," for values and peaks patterns.</p> <p>Note Leave a zero in the Size field to ignore a peak for the standard definition.</p>																								
c.	<p>Press Return to automatically move to the next size standard peak.</p>																								

To use the New command to define a new size standard: *(continued)*

Step	Action
8	<p>When you finish defining the peaks, save the size standard by choosing Save As from the File menu.</p> <p>Note You can also click the close box.</p>
9	<p>Type in a descriptive name for the standard and click Save.</p> <p>Note Run conditions are not stored in the Standard file. Use a name that clearly defines the standard for future use.</p> <p>This file is automatically saved in the Size Standards folder.</p> <p>For information on how to define the folder location, see page 5-41.</p>

Using the Analysis Control Window

To use the Analysis Control window to define a new size standard:

Step	Action					
1	<p>Open an existing project or create a new project.</p> <table border="1" data-bbox="467 808 991 927"> <thead> <tr> <th>For information on...</th> <th>See page</th> </tr> </thead> <tbody> <tr> <td>opening an existing project</td> <td rowspan="2">4-10</td> </tr> <tr> <td>creating a new project</td> </tr> </tbody> </table> <p>The Analysis Control window should appear. If it does not, choose Analysis Control (⌘ 1) from the Windows menu.</p>	For information on...	See page	opening an existing project	4-10	creating a new project
For information on...	See page					
opening an existing project	4-10					
creating a new project						
2	In the Analysis Control window, find the row that contains the sample for which you want to define the standard.					
3	<p>In that row, ⌘-click the dye color cell that represents your standard.</p> <p>A diamond symbol (◆) appears in the cell, identifying it as the standard.</p>					

To use the Analysis Control window to define a new size standard: *(continued)*

Step	Action								
4	<p>Click the arrow in the Parameters field of the same row, and select an option from the pop-up menu that appears.</p> <p>The pop-up menu contains the following options:</p> <table border="1" data-bbox="517 357 1229 820"> <thead> <tr> <th data-bbox="517 357 880 389">Item</th> <th data-bbox="880 357 1229 389">Description</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 389 880 495"><Analysis Parameters></td> <td data-bbox="880 389 1229 495">Applies the parameters that are stored as preferences in the software.</td> </tr> <tr> <td data-bbox="517 495 880 560"><filename></td> <td data-bbox="880 495 1229 560">Applies the settings specified in the data collection run file.</td> </tr> <tr> <td data-bbox="517 560 880 820">custom parameters that are listed at the bottom of the menu</td> <td data-bbox="880 560 1229 820"> <p>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p>For information on setting the Folder Locations preferences, see page 5-41.</p> </td> </tr> </tbody> </table>	Item	Description	<Analysis Parameters>	Applies the parameters that are stored as preferences in the software.	<filename>	Applies the settings specified in the data collection run file.	custom parameters that are listed at the bottom of the menu	<p>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p>For information on setting the Folder Locations preferences, see page 5-41.</p>
Item	Description								
<Analysis Parameters>	Applies the parameters that are stored as preferences in the software.								
<filename>	Applies the settings specified in the data collection run file.								
custom parameters that are listed at the bottom of the menu	<p>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences.</p> <p>For information on setting the Folder Locations preferences, see page 5-41.</p>								
5	<p>In the same row, click the arrow in the Size Standard field, and choose Define New from the pop-up menu.</p> <p>A window appears showing the electropherogram and a table of peaks for the dye color and sample you selected.</p> <p>You should be able to recognize the peak pattern of the standard in the electropherogram.</p>								
6	<p>Follow step 7 on page 5-33 to step 9 on page 5-34.</p> <p>The name of the standard appears in the following:</p> <ul style="list-style-type: none"> ◆ Standard pop-up menu in the Analysis Control window. ◆ Auto-Analysis settings dialog box. <p>You can specify in the Auto-Analysis Settings dialog box (or in the Data Collection software) that the defined standard be used for automatic analysis in future runs. See “Setting Up for Automatic Analysis” on page 4-4.</p>								

Using Size Standards

In This Section This section contains the following topics:

Topic	See page
Changing the Number of Peaks Detected	5-36
Editing the Size Standard Definition	5-36
Using the Open Command to Edit an Existing Standard	5-37
Using the Analysis Control Window to Edit an Existing Standard	5-38
Deleting an Existing Standard	5-39
Analyzing Samples Using the Same Standard	5-39
Selecting Separate Standards for Samples	5-40

Changing the Number of Peaks Detected

Use the Analysis Parameters dialog box to change the number of peaks detected in the Define New Standard window.

For information, see “Peak Detection Parameter Options” on page 5-20.

Editing the Size Standard Definition

The following procedure describes how to edit the size standard definition (<Collection Settings> standard) that is embedded in the Sample file. Double-clicking the size standard definition will not open the file.

To edit the size standard definition:


Step	Action
1	If the file is on your hard disk, choose Open from the file menu. The <Collection Settings> does not change.
2	Make any changes to the file.
3	Choose the edited file from the Standard pop-up menu to apply the changes.

Editing an Existing Standard

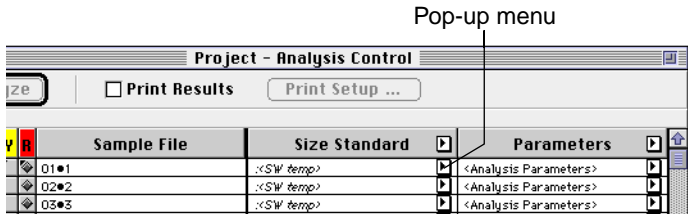
The two ways to edit a previously defined standard are by:

- ◆ Using the Open Command to Edit an Existing Standard.
- ◆ Using the Analysis Control Window to Edit an Existing Standard.

Using the Open Command to Edit an Existing Standard

Step	Action						
1	<p>Choose Open (⌘ O) from the File menu.</p> <p>The Open Existing dialog box appears.</p>						
	 <p>The screenshot shows a dialog box titled "Open Existing:". Inside, there are seven icons representing different file types: "Collection Gel" (spreadsheet), "Project" (project folder), "Sample" (graph), "Sample Sheet" (spreadsheet), "Analysis Parameters" (graph with a red dot), "Size Standard" (spreadsheet), and "Matrix" (grid). A "Cancel" button is located on the right side of the dialog box.</p>						
2	<p>Click the Size Standard icon.</p> <p>A directory dialog box appears.</p>						
3	<p>Select the standard file that you want to modify.</p>						
4	<p>Edit the peak size values by following step 7 to step 9 on page 5-34.</p>						
5	<p>Choose Save or Save As from the File menu.</p> <p>You can take the following action:</p> <table border="1" data-bbox="514 959 1239 1107"> <thead> <tr> <th>If you choose...</th> <th>Then the...</th> </tr> </thead> <tbody> <tr> <td>Save</td> <td>existing file is replaced.</td> </tr> <tr> <td>Save As</td> <td>changes are saved using a file name you specify.</td> </tr> </tbody> </table>	If you choose...	Then the...	Save	existing file is replaced.	Save As	changes are saved using a file name you specify.
If you choose...	Then the...						
Save	existing file is replaced.						
Save As	changes are saved using a file name you specify.						

Using the Analysis Control Window to Edit an Existing Standard

Step	Action												
1	If the Analysis Control window is not displayed, choose Analysis Control (⌘ 1) from the Windows menu.												
2	<p>Click the arrow in the Standard column for the sample that you want to change.</p> <p>A pop-up menu appears.</p>  <p>The screenshot shows the 'Project - Analysis Control' window. It has a search bar, a 'Print Results' checkbox, and a 'Print Setup ...' button. Below is a table with three columns: 'Sample File', 'Size Standard', and 'Parameters'. The 'Size Standard' column has a dropdown arrow for each row. A callout box labeled 'Pop-up menu' points to the dropdown arrow in the first row.</p> <table border="1"> <thead> <tr> <th>Sample File</th> <th>Size Standard</th> <th>Parameters</th> </tr> </thead> <tbody> <tr> <td>01#1</td> <td><:S\W temp></td> <td><Analysis Parameters></td> </tr> <tr> <td>02#2</td> <td><:S\W temp></td> <td><Analysis Parameters></td> </tr> <tr> <td>03#3</td> <td><:S\W temp></td> <td><Analysis Parameters></td> </tr> </tbody> </table>	Sample File	Size Standard	Parameters	01#1	<:S\W temp>	<Analysis Parameters>	02#2	<:S\W temp>	<Analysis Parameters>	03#3	<:S\W temp>	<Analysis Parameters>
Sample File	Size Standard	Parameters											
01#1	<:S\W temp>	<Analysis Parameters>											
02#2	<:S\W temp>	<Analysis Parameters>											
03#3	<:S\W temp>	<Analysis Parameters>											
3	Select a size standard from the pop-up menu.												
4	Edit the peak size values by following step 7 on page 5-33 to step 9 on page 5-34.												
5	<p>Choose Save or Save As from the File menu.</p> <p>You can take the following action:</p> <table border="1"> <thead> <tr> <th>If you choose...</th> <th>Then the...</th> </tr> </thead> <tbody> <tr> <td>Save</td> <td>existing file is replaced.</td> </tr> <tr> <td>Save As</td> <td>changes are saved using a file name you specify.</td> </tr> </tbody> </table>	If you choose...	Then the...	Save	existing file is replaced.	Save As	changes are saved using a file name you specify.						
If you choose...	Then the...												
Save	existing file is replaced.												
Save As	changes are saved using a file name you specify.												

Deleting an Existing Standard

The following procedure describes how to delete a user-defined standard from the GS Standards Folder so that it no longer appears in the Standard pop-up menus. The standard is permanently removed and you must re-define to use it again.

To delete an existing standard:

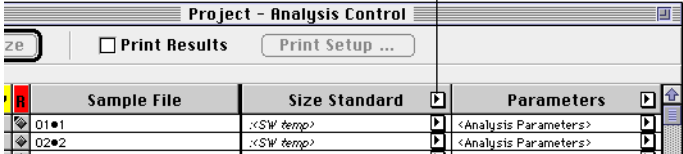
Step	Action
1	Switch to the Finder.
2	Locate and open the Current Standards folder.
3	Select the definition file that you want to delete.
4	Drag the file to the trash and empty the trash. Note You can also drag the standard to another folder for storage.

Analyzing Samples Using the Same Standard

To select the same standard for analysis of all samples:

Step	Action
1	If the Analysis Control window is not displayed, choose Analysis Control (⌘ 1) from the Windows menu.
2	Click the arrow in the Standard column heading and choose a standard file from the pop-up menu. Your menu choice applies to all fields in the column. Note Alternatively, you can choose a value from the pop-up menu, click the header to select the entire column, and choose Fill Down from the Edit menu.

Pop-up menu



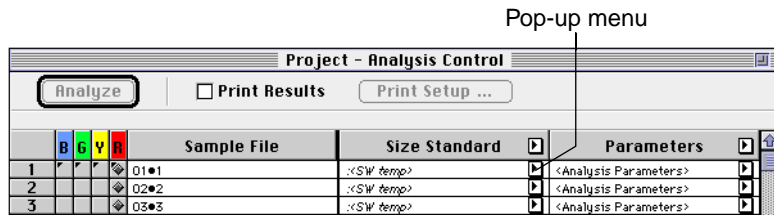
The screenshot shows the 'Project - Analysis Control' window. At the top, there are buttons for 'Print Results' and 'Print Setup ...'. Below is a table with three columns: 'Sample File', 'Size Standard', and 'Parameters'. The 'Size Standard' column has a pop-up menu open, showing a list of standard files. The table contains two rows of data, both with '01•1' and '02•2' in the 'Sample File' column and '<SW temp>' in the 'Size Standard' column. The 'Parameters' column contains '<Analysis Parameters>' for both rows.

Step	Action								
3	The pop-up menu contains the following options:								
	<table border="1"> <thead> <tr> <th>Item</th> <th>Description</th> </tr> </thead> <tbody> <tr> <td><None></td> <td>Apply no standard definition.</td> </tr> <tr> <td><filename></td> <td>Apply the size standard specified in the Data Collection software, which is embedded in the Sample file. For information on editing this file, see “Editing the Size Standard Definition” on page 5-36.</td> </tr> <tr> <td>custom standards that are listed at the bottom of the menu</td> <td>These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences. For information on setting the Folder Locations preferences, see “Defining Folder Locations” on page 5-41.</td> </tr> </tbody> </table>	Item	Description	<None>	Apply no standard definition.	<filename>	Apply the size standard specified in the Data Collection software, which is embedded in the Sample file. For information on editing this file, see “Editing the Size Standard Definition” on page 5-36.	custom standards that are listed at the bottom of the menu	These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences. For information on setting the Folder Locations preferences, see “Defining Folder Locations” on page 5-41.
	Item	Description							
	<None>	Apply no standard definition.							
<filename>	Apply the size standard specified in the Data Collection software, which is embedded in the Sample file. For information on editing this file, see “Editing the Size Standard Definition” on page 5-36.								
custom standards that are listed at the bottom of the menu	These are files that you defined and they are located in the parameters folder specified in the Folder Locations preferences. For information on setting the Folder Locations preferences, see “Defining Folder Locations” on page 5-41.								

Selecting Separate Standards for Samples

To apply separate size standards to selected samples, click the arrow in the Standard column for the sample that you want to change (see figure below) and a pop-up menu appears.

For information on using the pop-up menu, see step 3 above.



Defining Folder Locations

Introduction The GeneScan Analysis Software looks in the designated folders for the:

- ◆ Size Standard file.
- ◆ Analysis Parameter file.
- ◆ Matrix file.

When saving one of these files for the first time, the default folder locations for saving the files are those same designated folders.

Storing Matrix Files Store matrix files, that are intended for use by Data Collection software to assign to collection runs, in the ABI Folder. The ABI Folder is located in the System folder on the computer on which the Data Collection software is installed.

If Data Collection and Analysis are Performed on Different Computers

Make a copy of a matrix and store as follows. This is useful when data collection and analysis are performed on different computers.

Store a copy in the...	For use by the...
ABI Folder	Data Collection software.
GS Matrix Folder	GeneScan Analysis Software.

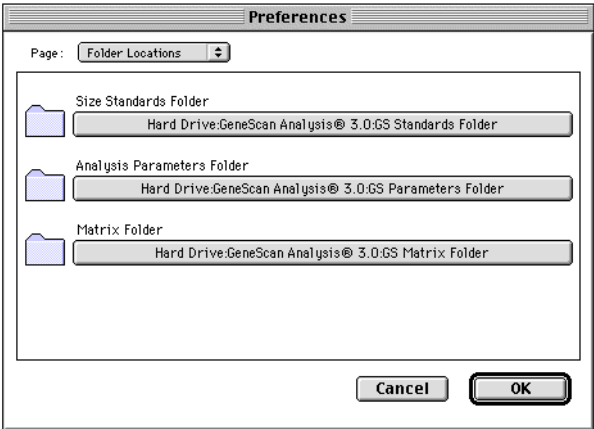
Note The ABI PRISM® instrument Data Collection software uses the files installed by the GeneScan Analysis Software in the ABI folder. When you run the analysis software, the program also creates several files (such as a Preference file) and an Analysis Log.

Specifying File Locations

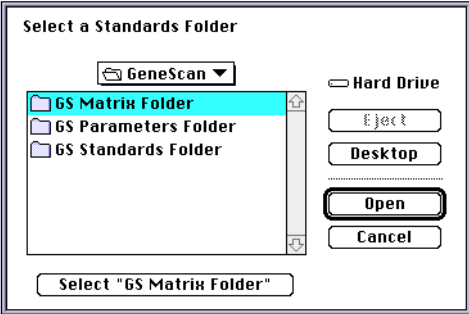
Use the Folder Location preference in the Settings menu to specify where the software saves size standards, analysis parameters, and matrices.

These preferences are saved as defaults for subsequent projects and Sample files.

To specify file locations:

Step	Action
1	<p>Choose Preferences from the Settings menu and Folder Locations from the submenu.</p> <p>The Preferences window appears (see below).</p> <p>Note It is advisable that you use the appropriate folders, that is, the GS Standards Folder, the GS Parameters Folder, and the GS Matrix Folder.</p> <p>Note If the Preference window is displayed, choose Folder Locations from the Page pop-up menu.</p>  <p>The screenshot shows a 'Preferences' dialog box with a 'Page:' dropdown menu set to 'Folder Locations'. Below the dropdown, there are three folder selection fields: 'Size Standards Folder' with the path 'Hard Drive:GeneScan Analysis@ 3.0:GS Standards Folder', 'Analysis Parameters Folder' with the path 'Hard Drive:GeneScan Analysis@ 3.0:GS Parameters Folder', and 'Matrix Folder' with the path 'Hard Drive:GeneScan Analysis@ 3.0:GS Matrix Folder'. At the bottom right of the dialog are 'Cancel' and 'OK' buttons.</p>

To specify file locations: *(continued)*

Step	Action						
2	<p>Click a button that shows the folder's path name. A dialog box appears.</p> 						
3	<p>Take the following action:</p> <ol style="list-style-type: none"> Find the new location. Highlight the target folder. Click the Select button at the bottom of the dialog box. 						
4	<p>Take the following action:</p> <table border="1" data-bbox="513 834 1239 1040"> <thead> <tr> <th data-bbox="513 834 876 873">If you have...</th> <th data-bbox="876 834 1239 873">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="513 873 876 971">other folder locations to designate</td> <td data-bbox="876 873 1239 971">choose another folder from the Folder Locations Preferences dialog box.</td> </tr> <tr> <td data-bbox="513 971 876 1040">no other folder locations to designate</td> <td data-bbox="876 971 1239 1040">click OK.</td> </tr> </tbody> </table>	If you have...	Then...	other folder locations to designate	choose another folder from the Folder Locations Preferences dialog box.	no other folder locations to designate	click OK.
If you have...	Then...						
other folder locations to designate	choose another folder from the Folder Locations Preferences dialog box.						
no other folder locations to designate	click OK.						

6

Making a Matrix File

Introduction

In This Chapter Topics in this chapter include the following:

Topics	See page
About Matrix Files	6-2
Process of Creating a New Matrix File	6-7
The Dye Matrix Standard Kits	6-8
Preparing Matrix Standards for the ABI PRISM 310	6-9
Preparing Matrix Standards for the ABI 373 and ABI PRISM 377	6-12
Loading and Running Dye Standards for the ABI PRISM 310	6-13
Loading and Running Dye Standards for the ABI PRISM 377	6-17
Loading and Running Matrix Samples on the ABI 373 (non XL)	6-20
Loading and Running Matrix Standards on the ABI 373 with XL Upgrade	6-24
Generating Matrix Sample Files for the ABI 373 and ABI PRISM 377	6-27
Choosing a Scan Range for the Matrix Calculation	6-29
Generating a New Matrix File	6-32
Saving and Naming the Matrix File	6-34
Assigning the Matrix File to Sample Files	6-35
Evaluating the Matrix File	6-37
Causes for Bad Matrix Files	6-38

About Matrix Files

Introduction There are three dye-labeling chemistries currently available to prepare nucleic acid samples for using the GeneScan® Analysis Software on ABI PRISM® instruments:

- ◆ Fluorescent NHS-Ester
- ◆ Fluorescent dNTP
- ◆ Fluorescent Phosphoramidite

Each chemistry has a set of dye labels that fluoresce at different wavelengths when excited by a laser.

During data collection on the...	The wavelengths are separated...
ABI PRISM® 310 or ABI PRISM® 377, 377XL, or 96-lane upgrade	by a spectrograph into a known spectral pattern across a detection system with the sequencer.
ABI™ 373 and the ABI™ 373 with XL upgrade	using a filter wheel.

Matrix File Definition Matrix files are mathematical matrices that correct for spectral overlap of fluorescent emission spectra data collected from ABI PRISM® instruments.

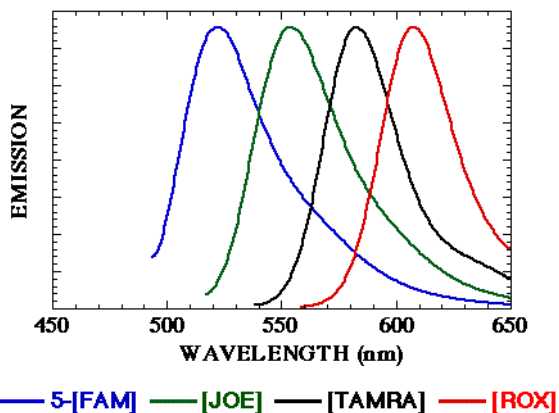
A matrix file allows you to account for spectral overlap when analyzing Sample files.

Multicomponent Definition This process of eliminating the bleed-through caused by spectral overlaps is called multicomponenting.

Applying a matrix file to raw data allows you to generate multicomponented data.

Why is a Matrix File Necessary

A matrix file is necessary because the four or five dyes used to label your fragments fluoresce at different wavelengths and may have spectral overlaps, as shown below.



When to Create a Matrix File

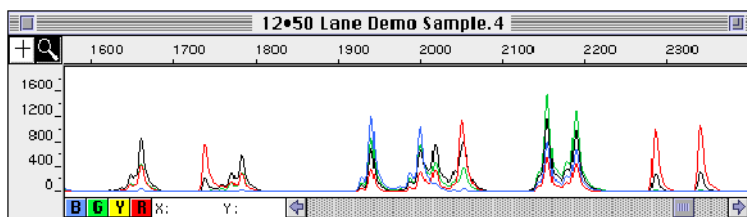
Create a matrix file for each dye set used from that particular instrument before analyzing fragment data.

You may have to create new matrix files for different gel compositions or unusual run conditions.

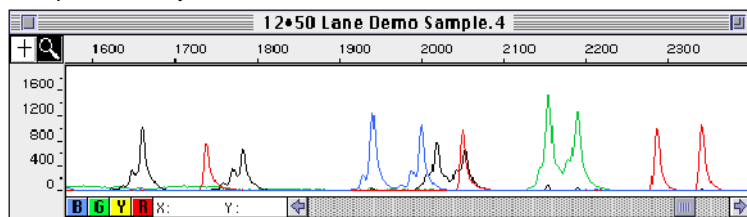
Sample Files Using Matrix File

The figure below shows an example of data analyzed with and without a matrix file.

You can see that peak data from a Sample file analyzed without a matrix file displays the expected peak, along with extra peaks in other dye colors, or bleed-through from other dye colors.



Sample file analyzed without a matrix file



Sample file analyzed with a matrix file

Applying a Matrix File

You can apply a matrix to a gel image or to the raw data within Sample files.

The matrix from a matrix file is installed within a gel file or Sample file automatically upon generation during or after a run, or manually from within the GeneScan Analysis Software.

If a matrix is installed to the...	See page
Gel image during image generation	2-24
Sample file's raw data during analysis	3-20

When to Assign a Matrix File

Before you can successfully analyze Sample files using the GeneScan Analysis Software, you must make a new matrix file or assign an existing one to a set of Sample files.

Limitations to Matrix Files

You can only assign a matrix file to Sample files generated on the same instrument, under the same electrophoresis, gel matrix and buffer conditions, and using the same dye set.

Note If you are using a fifth dye, then you need to create a new matrix file for that dye.

When to Create New a Matrix File

Create a new matrix file in the following conditions:

- ◆ For each dye set:
 - NHS-Esters
 - Phosphoramidite set
 - Fluorescent dNTPs
 - ◆ Whenever you change the dye set you use to label sample fragments, for example if you are using the fifth dye.
 - ◆ When you use gel materials or buffers with pH values that differ greatly from the pH value of the gel material or buffer on which the existing matrix files were generated.
 - ◆ When you use dyes other than those provided by PE Applied Biosystems.
 - ◆ When you run the same gel on a different instrument.
 - ◆ When you see multiple unexpected peaks of different colors under an expected peak.
 - ◆ When you recalibrate your CCD camera (ABI PRISM 310 and ABI PRISM 377) and the change is greater than 3 pixels from the original pixel position.
 - ◆ When you replace your filter wheel (ABI 373) or CCD camera (ABI PRISM 310 and ABI PRISM 377).
-
-

Considerations Before Making a Matrix File

The following table lists some of the considerations before making a matrix file:

Consideration	Comment
How much dye matrix standard to load?	With the ABI 373 or ABI PRISM 377, loading more than 3- μ L, produces too much signal. Any amount that results in a signal over 4,000 FU's is too strong.
Which lanes to load with the dye matrix standards?	For gel electrophoresis, load the matrix standards with an empty lane between each sample to avoid contamination of the individual dyes by residual material leaking adjacent samples.
After generating a gel image, for ABI 373 and ABI PRISM 377 instruments, check that the tracking of the gel file is adequate.	

Where to Store Matrix Files

Store matrix files intended for use by Data Collection software in the ABI folder. The ABI folder is located in the System folder on which the Data Collection software is installed.

If Data Collection and Analysis are Performed on Different Computers

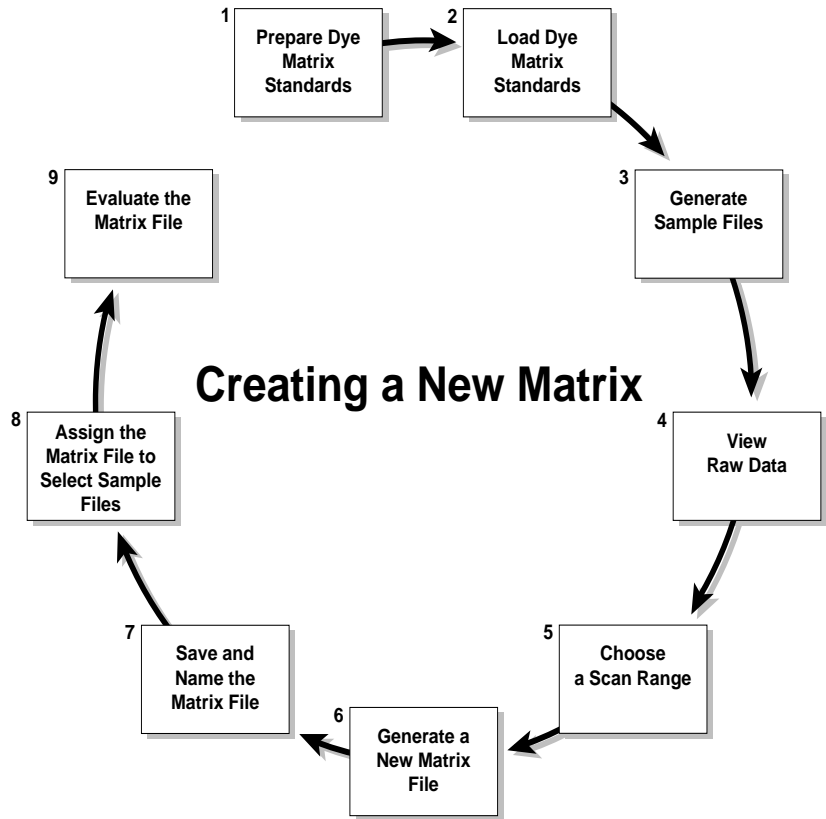
Make a copy of a matrix and store as follows. This is useful when data collection and analysis are performed on different computers.

Store a copy in the...	For use by the...
ABI Folder	Data Collection software.
GS Matrix Folder	GeneScan Analysis Software.

Note The ABI PRISM® instrument Data Collection software uses the files installed by the GeneScan Analysis Software in the ABI folder. When you run the analysis software, the program also creates several files (such as a Preference file) and an Analysis Log.

Process of Creating a New Matrix File

Process Diagram The following diagram shows the procedure for making a new matrix file:



Steps to Creating a New Matrix

The following table lists the steps to creating a new matrix file:

Step	Process	See page
1	Preparing Matrix Standards for the ABI PRISM 310.	6-9
2	Preparing Matrix Standards for the ABI 373 and ABI PRISM 377.	6-12
3	Loading and Running Dye Standards for the ABI PRISM 310.	6-13
4	Loading and Running Dye Standards for the ABI PRISM 377.	6-17
5	Loading and Running Matrix Samples on the ABI 373 (non XL).	6-20
6	Generating Matrix Sample Files for the ABI 373 and ABI PRISM 377.	6-27
7	Choosing a Scan Range for the Matrix Calculation.	6-29
8	Generating a New Matrix File.	6-32
9	Saving and Naming the Matrix File.	6-34
10	Assigning the Matrix File to Sample Files.	6-35
11	Evaluating the Matrix File.	6-37

The Dye Matrix Standard Kits

Table of Kits PE Applied Biosystems provides Dye Matrix Standard kits that are designed to test and define the multicomponent matrices for specific dye sets. Tubes containing matrix standards for each of the four different dye colors (blue, green, red, and yellow) are available in the each of these kits:

Kit	Part number
Dye Primer Matrix Standard Kit	401114
Fluorescent Amidite Matrix Standards	401456
Fluorescent dNTP Matrix Standards	402792

Preparing Matrix Standards for the ABI PRISM 310

About Formamide and Samples in Formamide

! WARNING ! CHEMICAL HAZARD. Formamide is a known teratogen. It can cause birth defects. Wash thoroughly after handling formamide. Wear appropriate protective eyewear, clothing, and gloves. Obtain a copy of the MSDS from the manufacturer. Wash thoroughly after handling formamide.

IMPORTANT Use the same matrix standards that you used for your GeneScan Analysis Software run.

The protocol uses formamide as a sample preparation reagent. Fresh formamide must be deionized and aliquotted into smaller volumes for storage. Each aliquot should be adequate for about one week's work.

Store aliquots of formamide at -20°C for up to three months. Formamide stored at 4°C is good for about one week.

At room temperature, samples in formamide are stable for a maximum of 48 hours.

Although not recommended on a routine basis, you can keep samples prepared in formamide frozen for no more than three days, with no detectable loss in resolution, before running on the ABI PRISM® 310 Genetic Analyzer.

Deionizing Formamide

To deionize formamide:

Step	Action	
1	Mix 50-mL of formamide and 5 g of ion-exchange resin (AG501 X8 from BioRad is recommended).	
2	Stir the mixture at room temperature for 30 minutes.	
3	Check the pH. It must be 7.0–9.0.	
4	If the pH...	Then...
	is in the proper range	filter the mixture through a 2-micron filter.
	is not in the proper range	a. Decant the formamide into a beaker with 5g of resin. b. Stir at room temperature for 30 minutes. c. Check the pH; it must be 7.0 – 9.0.

Step	Action
5	Make 500- μ L aliquots and store them at minus 20°C for up to 3 months.

Preparing the Formamide-Size Standard Mix

To ensure reproducibility of results for all samples, prepare the formamide-size standard mix using the 12:1 ratio of reagents stated in the procedure below.

IMPORTANT The formamide-size standard mix for the ABI PRISM 310 differs from the mix prepared for the ABI 373 and ABI PRISM 377, and do not use on other instruments.

! WARNING ! CHEMICAL HAZARD. Formamide is a known teratogen. It can cause birth defects. Wash thoroughly after handling formamide. Wear appropriate protective eyewear, clothing, and gloves. Obtain a copy of the MSDS from the manufacturer. Wash thoroughly after handling formamide.

Prepare the formamide-size standard mix as follows:

Step	Action
1	Mix in a sample vial: a. 0.5- μ L GeneScan-350 [TAMRA] Size Standard (for example). b. 12.0- μ L deionized formamide.
2	Label the vial.
3	Gently vortex the mixture for 3–5 seconds.
4	Spin down the mixture.
5	Store the mix at 2–6 °C until ready to use.

Preparing Matrix Standard Samples

About Matrix Standards

You must run matrix standards and create a matrix file the first time you use a new chemistry or change the run conditions.

Do not prepare matrix standards more than two hours in advance.

To ensure reproducibility of results for all samples, prepare the matrix standard mix using the 12:1 ratio of reagents stated in the procedure below.

IMPORTANT Do not add size standard mix to the matrix standard samples.

IMPORTANT The matrix standard mix for the ABI PRISM 310 differs from the mix prepared for the ABI 373 and ABI PRISM 377, and should not be used on other instruments.

For...	Prepare the...
the Fluorescent Genotyping Demonstration Kit	Fluorescent Amidite Matrix Standards (P/N 401546)
fluorescent dNTPs	Fluorescent dNTP Matrix Standards (P/N 402792).

Note Remember that the matrix standards run must always match the sample run chemistry and conditions.

Preparing Matrix Standard Samples

For each matrix standard:

Step	Action
1	Mix in a sample vial: a. 1.0- μ L of matrix standard. b. 12.0- μ L of de-ionized formamide.
2	Label each vial according to the dye.
3	Gently vortex the mixture for 3–5 seconds.
4	Spin down the mixture.
5	Store at 2–6 °C until ready to use.

Preparing Matrix Standards for the ABI 373 and ABI PRISM 377

Introduction Use the following procedure to prepare matrix standards for the following instruments:

ABI 373	ABI PRISM 377
ABI 373 with XL Upgrade	ABI PRISM® 377 with XL Upgrade

Procedure To prepare the matrix standards for loading:

Step	Action						
1	<p>Set up a run with the instrument conditions you want to duplicate.</p> <p>Run conditions depend on the following:</p> <ul style="list-style-type: none"> ◆ Instrument system you are using (ABI PRISM 377 or ABI 373). ◆ Kind of experiment that you are running. 						
2	<p>Select the four Dye Matrix Standards (one for each color) from the Dye Matrix Standard kit that you are using.</p> <table border="1" style="width: 100%;"> <thead> <tr> <th>If you are using...</th> <th>Then each dye that you use...</th> </tr> </thead> <tbody> <tr> <td>ABI PRISM 310 or ABI PRISM 377</td> <td>must be recognized as a different color by the single virtual filter.</td> </tr> <tr> <td>ABI 373</td> <td>must be recognized by a different filter on one filter wheel.</td> </tr> </tbody> </table>	If you are using...	Then each dye that you use...	ABI PRISM 310 or ABI PRISM 377	must be recognized as a different color by the single virtual filter.	ABI 373	must be recognized by a different filter on one filter wheel.
If you are using...	Then each dye that you use...						
ABI PRISM 310 or ABI PRISM 377	must be recognized as a different color by the single virtual filter.						
ABI 373	must be recognized by a different filter on one filter wheel.						
3	Depending on the instrument system, use the table below to prepare dye matrix standards for loading.						

Table of Matrix Standards The following table lists the matrix standards to prepare for each instrument:

For	ABI PRISM 377 and with XL Upgrade	ABI 373	ABI 373 with XL Upgrade
Denaturing gels	Combine 2.5- μ L of each standard with 1.5- μ L Loading buffer ^a .	Combine 2.5- μ L of each standard with 2.5- μ L Loading buffer.	Combine 2.5- μ L of standard with 2.5- μ L of deionized formamide.
Heating	Heat each sample at 95 °C for five minutes to denature before loading.		

a. Loading buffer is 1 part 25-mM EDTA/50-mg/ml Blue Dextran with 5 parts deionized formamide.

Loading and Running Dye Standards for the ABI PRISM 310

Introduction This section describes how to do the following tasks before running matrix standards on the ABI PRISM 310:

Topic	See page
Creating a GeneScan Sample Sheet	6-13
Creating a GeneScan Injection List	6-14
Starting the Data Collection Software	6-15

Note When loading the matrix standards on an instrument, note which colors you load in which autosampler positions.

Creating a GeneScan Sample Sheet

To create a GeneScan Sample Sheet:

Step	Action
1	Open the ABI PRISM 310 Data Collection software.
2	Choose New from the File menu.
3	Click the GeneScan Sample Sheet 48 or 96 tube icon. The GeneScan Sample Sheet appears.

Sample Name	Color	Std	Pres	Sample Info	Comments
A1 B (blue)	B	<input checked="" type="checkbox"/>	<input type="checkbox"/>		
	G	<input type="checkbox"/>	<input type="checkbox"/>		
	Y	<input type="checkbox"/>	<input type="checkbox"/>		
	R	<input type="checkbox"/>	<input type="checkbox"/>		
A3 G (green)	B	<input type="checkbox"/>	<input type="checkbox"/>		
	G	<input checked="" type="checkbox"/>	<input type="checkbox"/>		
	Y	<input type="checkbox"/>	<input type="checkbox"/>		
	R	<input type="checkbox"/>	<input type="checkbox"/>		
A5 Y (yellow)	B	<input type="checkbox"/>	<input type="checkbox"/>		
	G	<input type="checkbox"/>	<input type="checkbox"/>		
	Y	<input checked="" type="checkbox"/>	<input type="checkbox"/>		
	R	<input type="checkbox"/>	<input type="checkbox"/>		
A7 R (red)	B	<input type="checkbox"/>	<input type="checkbox"/>		
	G	<input type="checkbox"/>	<input type="checkbox"/>		
	Y	<input type="checkbox"/>	<input type="checkbox"/>		
	R	<input checked="" type="checkbox"/>	<input type="checkbox"/>		

4	Enter the appropriate colors (Blue, Green, Yellow, or Red) in the positions for each sample name.
---	---

To create a GeneScan Sample Sheet: *(continued)*

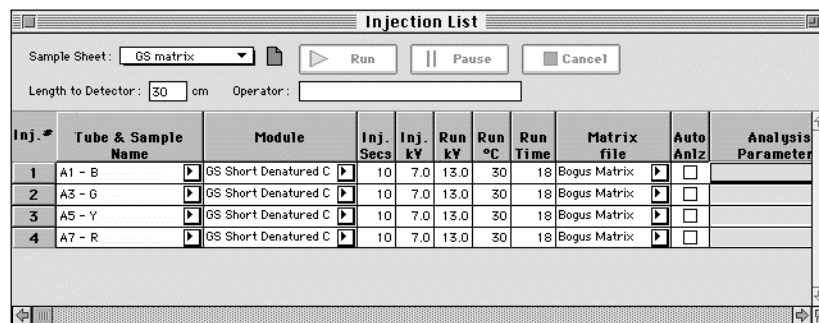
Step	Action
5	Enter any additional information about the sample in the Comments column.
6	Use the Save As command and save the Sample Sheet to the Sample Sheet folder.

Creating a GeneScan Injection List

The Sample Sheet is imported into the Injection List, which defines the sample names and the initial injection order.

To create a GeneScan Injection List:

Step	Action
1	Choose New from the File menu.
2	Click the GeneScan Injection List icon. The GeneScan Injection List appears.



3	Choose a Sample Sheet from the Sample Sheet pop-up menu.
4	Choose the appropriate module in the Module pop-up menu for lines 1 through 4, for example, A1, A3, A5, and A7.
5	Deselect Auto-Anlz (analysis).
6	Deselect Auto-Print.

Starting the Data Collection Software

To start the Data Collection software:

Step	Action						
1	Click the Run button. The following windows open:						
	<table border="1"> <thead> <tr> <th>Window</th> <th>Description</th> </tr> </thead> <tbody> <tr> <td>Raw Data window</td> <td>Shows the real-time chromatogram of the run.</td> </tr> <tr> <td>Log Window</td> <td>Shows the real-time written record of run events.</td> </tr> </tbody> </table>	Window	Description	Raw Data window	Shows the real-time chromatogram of the run.	Log Window	Shows the real-time written record of run events.
	Window	Description					
Raw Data window	Shows the real-time chromatogram of the run.						
Log Window	Shows the real-time written record of run events.						
2	Choose Status from the Window menu. The current run is in italics in the Injection List. You can monitor activities such as electrophoresis current, laser power, running time, and gel temperature.						

If the Run was Cancelled

The following table lists what happens to the Sample file if the run was cancelled:

If a run was cancelled and you are using...	Then the sample file is...
Data Collection software v 1.0.4 Note Version 1.0.4 is a free upgrade. To obtain the upgrade, either access the ABI PRISM 310 Genetic Analyzer website at http://www.perkin-elmer.com/ga/310 or call technical support (see page 1-20). For more information, refer to <i>ABI PRISM 310 Data Collection Software Version 1.0.4 User Bulletin</i> (P/N 4305798).	saved if you skip to the next sample or cancel a run.

Run Time Run time is approximately 30 minutes for the GS STR POP-4 module, so the total run time will be about 120 minutes.

Assigning Matrix File For information on Assigning the Matrix File to Sample Files, see page 6-35.

Loading and Running Dye Standards for the ABI PRISM 377

Introduction This section describes how to do the following tasks using the ABI PRISM 377 and the ABI PRISM 377 with XL Upgrade:

Topic	See page
Creating a GeneScan Sample Sheet	6-17
Loading Matrix Standards	6-18
Running the Matrix Standards	6-19

Note When loading the matrix standards on an instrument, note which colors you load in which lanes for gel based systems.

Creating a GeneScan Sample Sheet

The GeneScan Sample Sheet assigns sample and dye information to their appropriate lane.

To create a GeneScan Sample Sheet:

Step	Action
1	Open the ABI PRISM 377 Data Collection software. Define or verify the data collection preferences.
2	Choose New from the File menu.
3	Click the GeneScan Sample Sheet icon. The GeneScan Sample Sheet appears.

Sample Sheet "GS matrix"						
GeneScan™ Sample Sheet						
#	Sample Name	Color	Std	Pres	Sample Info	Comments
A1	B (blue)	B		<input checked="" type="checkbox"/>		
		G		<input type="checkbox"/>		
		Y		<input type="checkbox"/>		
		R		<input type="checkbox"/>		
A3	G (green)	B		<input type="checkbox"/>		
		G		<input checked="" type="checkbox"/>		
		Y		<input type="checkbox"/>		
		R		<input type="checkbox"/>		
A5	Y (yellow)	B		<input type="checkbox"/>		
		G		<input type="checkbox"/>		
		Y		<input checked="" type="checkbox"/>		
		R		<input type="checkbox"/>		
A7	R (red)	B		<input type="checkbox"/>		
		G		<input type="checkbox"/>		
		Y		<input type="checkbox"/>		
		R		<input checked="" type="checkbox"/>		

To create a GeneScan Sample Sheet: *(continued)*

Step	Action
4	Enter the individual colors in the appropriate lanes where the matrix standards are loaded. Note It is important to fill out the Sample Sheet completely.
5	Enter any additional information about the sample in the Comments column.
6	Use the Save As command and save the Sample Sheet to the Sample Sheet folder.

Loading Matrix Standards

To load matrix standards:

Step	Action									
1	For denaturing gels load: <ul style="list-style-type: none"> ◆ 0.5-2-μL of matrix standard per lane. ◆ 8 lanes with different colors, leaving an empty lane between each lane of matrix standard. 									
2	Complete the information in the data collection Run sheet, making sure to choose the appropriate PreRun and Run modules. Take the following action: <table border="1" style="margin-left: 20px;"> <thead> <tr> <th>For this matrix...</th> <th>Choose modules that use...</th> <th>Module file</th> </tr> </thead> <tbody> <tr> <td>Dye Primer matrix</td> <td>Virtual Filter A</td> <td>GS 36A...</td> </tr> <tr> <td>Fluorescent Amidite</td> <td>Virtual Filter C</td> <td>GS36C...</td> </tr> </tbody> </table>	For this matrix...	Choose modules that use...	Module file	Dye Primer matrix	Virtual Filter A	GS 36A...	Fluorescent Amidite	Virtual Filter C	GS36C...
For this matrix...	Choose modules that use...	Module file								
Dye Primer matrix	Virtual Filter A	GS 36A...								
Fluorescent Amidite	Virtual Filter C	GS36C...								
3	Electrophorese samples according to conditions specified in your instrument manual.									

Running the Matrix Standards

Run the matrix standards under the precise conditions you want to generate a matrix file.

To run the matrix standards:

Step	Action
1	Complete the information in the data collection Run sheet. Refer to the figure below to choose the appropriate Prerun and Run modules.

Lane	Sample Number	Sample Name	Sample File Name	Matrix	Auto Analyze	Analysis Parameter
1				<none>	<input type="checkbox"/>	<none>
2				<none>	<input type="checkbox"/>	<none>
3				<none>	<input type="checkbox"/>	<none>
4				<none>	<input type="checkbox"/>	<none>
5				<none>	<input type="checkbox"/>	<none>

2	Take the following action:									
	<table border="1"> <thead> <tr> <th>For this matrix...</th> <th>Choose modules that use...</th> <th>Module file</th> </tr> </thead> <tbody> <tr> <td>Dye Primer or dNTP matrix</td> <td>Virtual Filter A</td> <td>GS 36A...</td> </tr> <tr> <td>GS Amidite matrix</td> <td>Virtual Filter C</td> <td>GS36C...</td> </tr> </tbody> </table>	For this matrix...	Choose modules that use...	Module file	Dye Primer or dNTP matrix	Virtual Filter A	GS 36A...	GS Amidite matrix	Virtual Filter C	GS36C...
For this matrix...	Choose modules that use...	Module file								
Dye Primer or dNTP matrix	Virtual Filter A	GS 36A...								
GS Amidite matrix	Virtual Filter C	GS36C...								
3	Start the electrophoresis run according to the conditions specified in your instruction manual.									
4	Go to "Generating Matrix Sample Files for the ABI 373 and ABI PRISM 377" on page 6-27.									

Loading and Running Matrix Samples on the ABI 373 (non XL)

Introduction The ABI 373 uses the 672 Data Collection software. For analysis of the data after completing the matrix run, transfer the data collection file to a Power Macintosh® computer.

This section describes how to do the following tasks:

Topic	See page
Loading the Matrix Standards	6-20
Completing the Sample Sheet in the 672 Software	6-21
Running the Gel	6-22
Completing the GeneScan Sample Sheet	6-23

Note See Appendix A for converting ABI 373 data collection data into gel files.

Loading the Matrix Standards

To load matrix standards:

Step	Action						
1	For denaturing and native gels load: <ul style="list-style-type: none"> ◆ 1-5-µL of matrix standard per lane (depending on comb configuration, see instrument user's manual). ◆ 8 lanes with different colors, leaving an empty lane between each lane of matrix standard. 						
2	Select the appropriate filter set on the ABI 373 sequencer. <table border="1" style="margin-left: 20px;"> <thead> <tr> <th>For this matrix...</th> <th>Choose...</th> </tr> </thead> <tbody> <tr> <td>Dye Primer matrix</td> <td>Filter set A.</td> </tr> <tr> <td>Fluorescent Amidite matrix</td> <td>Filter set B.</td> </tr> </tbody> </table>	For this matrix...	Choose...	Dye Primer matrix	Filter set A.	Fluorescent Amidite matrix	Filter set B.
For this matrix...	Choose...						
Dye Primer matrix	Filter set A.						
Fluorescent Amidite matrix	Filter set B.						
3	Electrophorese samples according to conditions specified in your instrument manual.						

Completing the Sample Sheet in the 672 Software

Introduction

After opening the Collection file, complete the Sample Sheet embedded and track the lanes. This ensures that when the Sample files are generated they contain the correct information.

The Sample Sheet that is part of the Collection file is blank when it is opened. This is because the GeneScan® 672 Software does not automatically embed a copy of the data collection Sample Sheet in the Collection file.

Procedure

IMPORTANT When tracking and extracting the Collection file, ensure the Sample Sheet associated with the gel file has the checkbox labeled “Used” selected for each sample you want to extract. The GeneScan Analysis Software only extracts the samples with that checkbox selected.

To complete the Sample Sheet:

Step	Action								
1	Open the Sample Sheet by either: <ul style="list-style-type: none"> ◆ Clicking the Sample Sheet button in the upper-left corner of the Gel Window, or by ◆ Choosing Gel Sample Sheet from the Gel menu. 								
2	Enter the following information: <ul style="list-style-type: none"> ◆ Sample file name. ◆ Sample name for each sample. ◆ Sample info. ◆ Comment for each dye/sample. For more information, see “Displaying the Sample Sheet” on page 2-29.								
3	Select checkboxes, as applicable. <table border="1" style="width: 100%; margin-top: 10px;"> <thead> <tr> <th>To...</th> <th>Then select the...</th> </tr> </thead> <tbody> <tr> <td>indicate that the lane contains data</td> <td>Used checkbox.</td> </tr> <tr> <td>automatically analyze Sample files when generating them</td> <td>"A" checkbox.</td> </tr> <tr> <td>automatically print Sample files when generating them</td> <td>"P" checkbox.</td> </tr> </tbody> </table>	To...	Then select the...	indicate that the lane contains data	Used checkbox.	automatically analyze Sample files when generating them	"A" checkbox.	automatically print Sample files when generating them	"P" checkbox.
To...	Then select the...								
indicate that the lane contains data	Used checkbox.								
automatically analyze Sample files when generating them	"A" checkbox.								
automatically print Sample files when generating them	"P" checkbox.								

To complete the Sample Sheet: *(continued)*

Step	Action
4	Identify the dye standard for each sample by clicking in the appropriate dye cells under the standard column. Note Alternately, you can create a stand-alone Sample sheet and attach it to the gel file using the Install New Sample Sheet command from the Gel menu (see "Installing a New Sample Sheet" on page 2-36).
5	Choose Save (⌘-S) or Save As from the File menu. If you click the close box before saving, a dialog box appears with the message whether or not to save the entries.

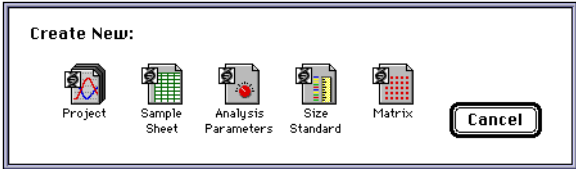
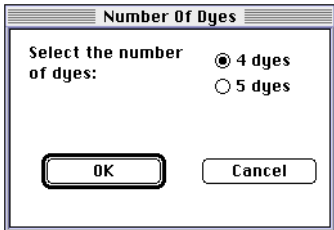
Running the Gel

To run the gel:

Step	Action						
1	Open the Data Collection software.						
2	Set the run time and Data Collection File name.						
3	On the ABI 373 instrument, choose the appropriate filter set for the conditions you wish to duplicate: <table border="1" data-bbox="481 816 1056 937" style="margin-left: 40px;"> <thead> <tr> <th>For this matrix...</th> <th>Choose...</th> </tr> </thead> <tbody> <tr> <td>Dye Primer or dNTP matrix</td> <td>Filter Set A.</td> </tr> <tr> <td>Fluorescent Amidite matrix</td> <td>Filter Set B.</td> </tr> </tbody> </table>	For this matrix...	Choose...	Dye Primer or dNTP matrix	Filter Set A.	Fluorescent Amidite matrix	Filter Set B.
For this matrix...	Choose...						
Dye Primer or dNTP matrix	Filter Set A.						
Fluorescent Amidite matrix	Filter Set B.						
4	Run the matrix standards.						
5	Start the electrophoresis run according to the conditions specified in your instruction manual.						

Completing the GeneScan Sample Sheet

To complete the Sample Sheet in the GeneScan Analysis Software:

Step	Action								
1	Open the GeneScan Analysis Software program.								
2	Choose New from the File menu. The Create New dialog box appears. 								
3	Click the Sample Sheet icon. The following dialog box appears: 								
4	Click the radio button that corresponds to the number of dyes and click OK. An untitled Sample Sheet appears.								
5	Fill out the Sample Sheet for each of the colors.								
6	Set the preprocessing parameters as follows: <table border="1" data-bbox="514 1065 1239 1284"> <thead> <tr> <th>Do...</th> <th>And...</th> </tr> </thead> <tbody> <tr> <td>deselect Multicomponent</td> <td></td> </tr> <tr> <td>select Use Sample Sheet</td> <td>choose the appropriate Sample Sheet.</td> </tr> <tr> <td>select Auto Lane Track Baseline Data</td> <td>leave the Process amount as the default value.</td> </tr> </tbody> </table>	Do...	And...	deselect Multicomponent		select Use Sample Sheet	choose the appropriate Sample Sheet.	select Auto Lane Track Baseline Data	leave the Process amount as the default value.
Do...	And...								
deselect Multicomponent									
select Use Sample Sheet	choose the appropriate Sample Sheet.								
select Auto Lane Track Baseline Data	leave the Process amount as the default value.								
7	Choose Install New Sample Sheet from the Gel menu.								
8	Go to "Generating Matrix Sample Files for the ABI 373 and ABI PRISM 377" on page 6-27.								

Loading and Running Matrix Standards on the ABI 373 with XL Upgrade

Introduction The ABI 373 with XL upgrade updates the Data Collection software of the ABI 373 to make it compatible with Power Macintosh® computers, and allows you to open gel files directly using the GeneScan Analysis Software in much the same way the ABI PRISM 377 does.

This section describes how to do the following tasks:

Topic	See page
Completing the GeneScan Sample Sheet	6-24
Loading the Matrix Standards	6-25
Running the Matrix Standard	6-26

Note The procedures in Appendix A are not necessary since the ABI 373 software operates on a Power Macintosh computer.

Completing the GeneScan Sample Sheet

The GeneScan Sample Sheet assigns the sample and dye information to their appropriate lanes.

To create a GeneScan Sample Sheet:

Step	Action
1	Open the Data Collection software.
2	Choose New from the File menu.
3	Click the GeneScan Sample Sheet icon. The GeneScan Sample Sheet appears.

To create a GeneScan Sample Sheet: *(continued)*

Step	Action
4	Enter the individual colors in the appropriate lanes where the matrix standards are loaded.
5	Enter any additional information about the sample in the Comments column.
6	Use the Save As command and save the Sample Sheet to the Sample Sheet folder.

Loading the Matrix Standards

To load matrix standards:

Step	Action						
1	<p>For denaturing and native gels load:</p> <ul style="list-style-type: none"> ♦ 1-5-μL of matrix standard per lane (depending on comb configuration, see instrument user's manual). ♦ 8 lanes with different colors, leaving an empty lane between each lane of matrix standard. 						
2	<p>Select the appropriate filter set on the ABI 373 sequencer.</p> <table border="1" style="width: 100%; margin-top: 10px;"> <thead> <tr> <th>For this matrix...</th> <th>Choose...</th> </tr> </thead> <tbody> <tr> <td>Dye Primer matrix</td> <td>Filter set A.</td> </tr> <tr> <td>Fluorescent Amidite matrix</td> <td>Filter set B.</td> </tr> </tbody> </table>	For this matrix...	Choose...	Dye Primer matrix	Filter set A.	Fluorescent Amidite matrix	Filter set B.
For this matrix...	Choose...						
Dye Primer matrix	Filter set A.						
Fluorescent Amidite matrix	Filter set B.						

Step	Action
3	Electrophorese samples according to conditions specified in your instrument manual.

Running the Matrix Standard

Run the matrix standards under the precise conditions you want to generate a matrix file.

To run the matrix standards:

Step	Action
1	Complete the information in the Run window. Note Make sure to choose the appropriate PreRun and Run modules.

2	Choose the appropriate filter set for the conditions you wish to duplicate. <table border="1" data-bbox="467 1214 1193 1333"> <thead> <tr> <th>For this matrix...</th> <th>Choose...</th> </tr> </thead> <tbody> <tr> <td>Dye Primer or dNTP matrix</td> <td>Filter Set A.</td> </tr> <tr> <td>Fluorescent Amidite matrix</td> <td>Filter Set B.</td> </tr> </tbody> </table>	For this matrix...	Choose...	Dye Primer or dNTP matrix	Filter Set A.	Fluorescent Amidite matrix	Filter Set B.
For this matrix...	Choose...						
Dye Primer or dNTP matrix	Filter Set A.						
Fluorescent Amidite matrix	Filter Set B.						
3	Start the electrophoresis run according to the conditions specified in your instruction manual.						

Generating Matrix Sample Files for the ABI 373 and ABI PRISM 377

Who Should Use This Step

Who should use this step	Who shouldn't use this step
<p>This step is only necessary if you are using an:</p> <ul style="list-style-type: none"> ◆ ABI 373 ◆ ABI PRISM 377 ◆ XL-upgraded instrument ◆ 96-lane upgrade 	<p>The ABI PRISM 310 Genetic Analyzer automatically processes collection data and generates Sample files when the run completes.</p> <p>If you ran your Dye Matrix Standards on the ABI PRISM 310, go to "Choosing a Scan Range for the Matrix Calculation" on page 6-29.</p>

Verify Tracking

Before generating Sample files, verify that the lanes were set and tracked correctly. The following table lists where the data appears for each instrument once you have successfully completed your run of Dye Matrix Standards.

On this instrument...	The data appears...	Next step...
ABI PRISM 377	in a Gel File window.	Verify that the lanes were set and tracked correctly.
ABI 373	in a Data Collection file. Convert the Data Collection file to generate the gel image (see Appendix A).	

Generating Sample Files

To generate Sample files:

Step	Action
1	<p>Click the appropriate Lane Indicators at the top of the Gel window to make sure that the tracker line is optimally aligned over each band in all lanes.</p> <p>If any of the lanes are not properly tracked, use the tracker line editing tools to align the tracker lines in each lane. See "Working with Tracker Lines" on page 2-53.</p>

To generate Sample files: *(continued)*

Step	Action
2	If you change any of the lane assignments in any way, save the changes.
3	From the Gel menu, choose Track & Extract Lanes. Note If the gel is already tracked, choose Extract Lanes from the Gel menu. The project Analysis Control window appears containing each of the Sample files.

Choosing a Scan Range for the Matrix Calculation

Introduction Depending on how well your Matrix Standards run, it may be necessary for you to choose a specific range of data points to be considered for your matrix calculation.

In order to choose appropriate values for the Scan range, you must first view the Sample file raw data from each of the four matrix standard files, so you can decide where to choose the start and stop points for the scan range.

Viewing the Raw Data When you have multiple Sample files, raw data can be accessed more easily through a project's Analysis Control window. Raw data can provide useful information about the Sample files you have created.

Note You can view Sample files without opening a project. However, this procedure is easier if you use a project to organize the Sample files (see "Using a Project to Manage Sample Files" on page 4-8).

To view raw data:

Step	Action														
1	Use the following steps to create a project for the Dye Matrix Standards:														
	<table border="1"><thead><tr><th>Step</th><th>Action</th></tr></thead><tbody><tr><td>a.</td><td>Choose New from the File menu. The Create New dialog box appears.</td></tr><tr><td>b.</td><td>Click the Project icon. An untitled Analysis Control window opens.</td></tr><tr><td>c.</td><td>Choose Add Sample Files from the Project menu.</td></tr><tr><td>d.</td><td>Find and open your matrix run folder.</td></tr><tr><td>e.</td><td>Select the four Sample files representing the blue, green, yellow, and red dye-labeled runs, and then click Add.</td></tr><tr><td>f.</td><td>Click Done after the Sample files are transferred.</td></tr></tbody></table>	Step	Action	a.	Choose New from the File menu. The Create New dialog box appears.	b.	Click the Project icon. An untitled Analysis Control window opens.	c.	Choose Add Sample Files from the Project menu.	d.	Find and open your matrix run folder.	e.	Select the four Sample files representing the blue, green, yellow, and red dye-labeled runs, and then click Add.	f.	Click Done after the Sample files are transferred.
	Step	Action													
	a.	Choose New from the File menu. The Create New dialog box appears.													
	b.	Click the Project icon. An untitled Analysis Control window opens.													
	c.	Choose Add Sample Files from the Project menu.													
	d.	Find and open your matrix run folder.													
	e.	Select the four Sample files representing the blue, green, yellow, and red dye-labeled runs, and then click Add.													
f.	Click Done after the Sample files are transferred.														
For more information, see "Creating a New Project" on page 4-10.															

To view raw data: *(continued)*

Step	Action
2	From the Analysis Control window, select the four matrix standard Sample files by clicking on the first Sample file, holding down the mouse button, and releasing on the last Sample file.
3	<p>From the Project menu, choose Raw Data (⌘ R).</p> <p>Electropherograms displaying raw data from the four matrix standard Sample files appear.</p> <p>Note For the ABI PRISM 377 or ABI 373, you can also view raw data from the gel display by selecting one of the four lanes containing Dye Matrix Standard and looking at the Slice View to the left of the gel image.</p> <p>For more information about viewing the gel image, see “What to Review in the Gel Image” on page 2-27.</p>

What to Look For in the Raw Data Display

In the raw data display of the Sample files verify the following:

- ◆ Data peaks are present in all four of the matrix standards.
- ◆ There are no anomalies.
- ◆ The baseline is stable.
- ◆ Peaks should be on-scale—no more than 4,000 relative fluorescent units—and the peaks of the dye of interest should have a value of at least 200.

If peak data does not show these characteristics, see “Causes for Bad Matrix Files” on page 6-38, for possible interpretations of your peak data.

Choosing a Scan Range

To choose a scan range:

Step	Action
1	<p>Move the cursor well away from the primer peak, in a region at the beginning of the run and in a flat part of the baseline, and record the scan numbers.</p> <p>Note When choosing the start point do not include primer peaks in the scan range (see “Eliminating Primer Peaks” on page 6-31). Also, the region for both the start and stop points should be flat points at the baseline.</p>

Step	Action
2	Record the data point values for both the start and stop points of the scan range. You will need to enter those values in the next step when generating the new matrix file (see page 6-32).
3	Close the raw data boxes and the project by clicking in the upper left-hand corner of the window. Note Holding down the Option key while clicking in the upper left-hand corner of the window will close the windows simultaneously.

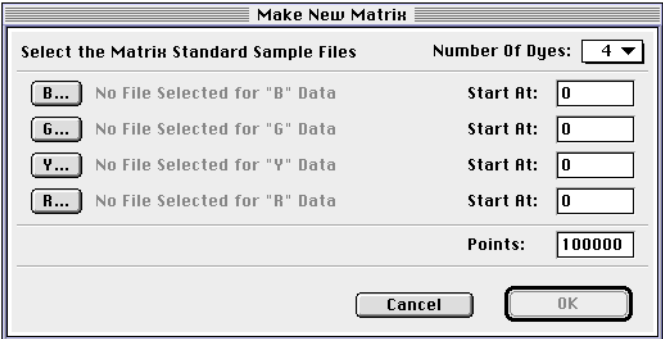
Eliminating Primer Peaks

Both the primer peaks and the data peaks are displayed when viewing the raw data of your matrix standards. Anytime you run dye-labeled samples on a gel (ABI 373 and ABI PRISM 377), or capillary (ABI PRISM 310), you always have excess dye-labeled primer in the reaction. The primer peak displays as the first peak, usually off-scale because it is in molar excess.

Eliminate the primer peak when making a matrix by choosing the start point after the primer peak in a flat area with a stable baseline.

Generating a New Matrix File

Procedure To generate a new matrix file:

Step	Action
1	Choose New from the File menu. The Create New dialog box appears.
2	Click the Matrix icon. The Make New Matrix dialog box appears. 
3	Choose the number of dyes from the Number of Dyes pop-up menu. If 5 dyes are selected, a button is added to the bottom of the list.
4	The B, G, Y, and R buttons represent dye colors. <ul style="list-style-type: none"> a. Click a button to display a pop-up menu. b. Use the pop-up menu to access a Sample file to link to each of the four dye-labeled primers. c. Choose the Sample file that represents the dye color for that button.
5	Enter the start point that you determined when choosing a scan range in the Start at field. See "Choosing a Scan Range" on page 6-30.
6	Enter the total number of data points to include to calculate the matrix in the Points field. <p>Note You must have at least five peaks to make a matrix.</p> <p>In most cases, leave the default value, unless you must exclude a portion of your data because of artifacts or bleed-through.</p>

Step	Action
7	Click OK. This generates a new matrix file.

Matrix File Example The following is an example of the Matrix Values window that appears showing the values used to calculate the overlap correction.

The screenshot shows a window titled "Run Matrix" with a sub-header "Reactions". Below this is a 4x4 grid of input fields. The columns are labeled B, G, Y, and R. The rows are labeled B, G, Y, and R. Each cell contains a numerical value representing the matrix value for that reaction combination.

	B	G	Y	R
B	1.0000	0.2304	0.0073	0.0019
G	0.5916	1.0000	0.2640	0.0028
Y	0.3173	0.5916	1.0000	0.1445
R	0.1211	0.2663	0.5000	1.0000

For each dye, the value where the dye fluorescence is read by the appropriate filter is 1.000. The adjacent colors show the amount of overlap for which the system must compensate. The adjacent values, in most cases, should be less than 1.000, but equal to or greater than 0.0000.

Note In some Filter Set C matrices, the matrix value for green in the first column can vary from about 0.8 to 1.2.

Saving and Naming the Matrix File

Introduction The matrix file is instrument-specific. You cannot apply a matrix file you made on the ABI 373 to data you collected on a ABI PRISM 377 or ABI PRISM 310, nor can you apply a matrix file made on a ABI 373 to a Sample or gel file made on another ABI 373. Also, you cannot apply matrix files created on one instrument to other instruments of the same model.

Naming Considerations When naming a matrix file, consider including the following information in the name:

Item	For example
Instrument type	ABI 373, ABI PRISM 377, or ABI PRISM 310.
Filter set used	A, B, or C.
Gel conditions	native or denaturing.

Saving the Matrix File To save the matrix file:

Step	Action
1	If a matrix folder does not already exist inside your GeneScan applications folder, create one.
2	Choose Save from the File menu. A dialog box appears.
3	Enter a descriptive name for the new matrix file and click Save.

Where to Store the Matrix File Store matrix files that are intended for use by data collection to assign to collection runs in the ABI folder. The ABI folder is located in the System folder of the computer on which Data Collection software is installed.

When to Make Copies of a Matrix You can make a copy of a matrix and store one in the ABI Folder for use by the Data Collection software and the other in the GS Matrix Folder for use by the GeneScan Analysis Software. This is useful when data collection and analysis are performed on different computers.

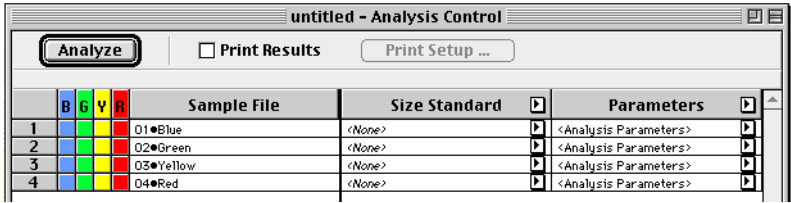
Assigning the Matrix File to Sample Files

Introduction After generating the new matrix file, assign it to all the Sample files that you want to analyze.

IMPORTANT After assigning your matrix file to Sample files, see “Evaluating the Matrix File” on page 6-37.

Procedure

To assign a matrix file to Sample files:

Step	Action
Assigning a matrix file to Sample files:	
1	From the project that contains your matrix standard Sample files, open the Analysis Control window.
	
2	Select the four matrix standard Sample files.
3	Choose Assign New Matrix from the Project menu.
4	Select the matrix file you just created.
5	Select numbers 1, 2, 3, and 4 on the left side of the window to highlight the colors for each row.
6	Choose Set Analysis Parameters from the Settings menu.
7	Enter the appropriate range for the Analysis Range and click OK to return to the Analysis Control window.
8	Click Analyze.
Reviewing the results:	
1	Choose Results Control from the Windows menu. The Results Control window appears (see below).

To assign a matrix file to Sample files: *(continued)*

Step	Action
2	Select 4 from the # of Panels pull-down menu (see above).
3	Click 1 under Dye/Samples (see above).
4	Click 1 on the Sample Files side of the Results control window.
5	For matrices 2, 3, and 4,click Analyze and repeat step 1 and step 2.

Evaluating the Matrix File

Introduction After creating a new matrix file and assigning it to select Sample files, the next step is evaluating the quality of the matrix file. The quality of the matrix files has a direct impact on the quality of the results data.

Procedure To evaluate the matrix file:

Step	Action						
1	Analyze the Sample files used to make the matrix.						
2	Display Results data for all four Dye Matrix Standard Sample files on one screen, showing only electropherogram data.						
3	For each displayed Sample file: <table border="1"><thead><tr><th>You should see...</th><th>If not...</th></tr></thead><tbody><tr><td>that the only visible peaks represent the color of the Dye Matrix Standard run in that lane, or for that injection (ABI PRISM 310).</td><td rowspan="3">you probably have a bad matrix file. For instructions on how to identify and correct problems with bad matrix files see, "Causes for Bad Matrix Files" on page 6-38.</td></tr><tr><td>all other lines should be relatively flat. This indicates that the matrix properly compensated for the spectral overlap. For example, for the blue matrix standard Sample file, you should only see blue.</td></tr><tr><td>sharp, well-defined, singularly colored peak data.</td></tr></tbody></table>	You should see...	If not...	that the only visible peaks represent the color of the Dye Matrix Standard run in that lane, or for that injection (ABI PRISM 310).	you probably have a bad matrix file. For instructions on how to identify and correct problems with bad matrix files see, "Causes for Bad Matrix Files" on page 6-38.	all other lines should be relatively flat. This indicates that the matrix properly compensated for the spectral overlap. For example, for the blue matrix standard Sample file, you should only see blue.	sharp, well-defined, singularly colored peak data.
You should see...	If not...						
that the only visible peaks represent the color of the Dye Matrix Standard run in that lane, or for that injection (ABI PRISM 310).	you probably have a bad matrix file. For instructions on how to identify and correct problems with bad matrix files see, "Causes for Bad Matrix Files" on page 6-38.						
all other lines should be relatively flat. This indicates that the matrix properly compensated for the spectral overlap. For example, for the blue matrix standard Sample file, you should only see blue.							
sharp, well-defined, singularly colored peak data.							

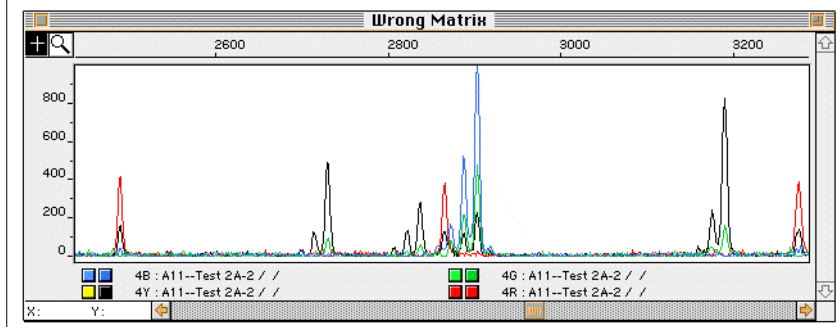
Causes for Bad Matrix Files

If an Error Message Appears If an error message appears, the following table lists two possible causes:

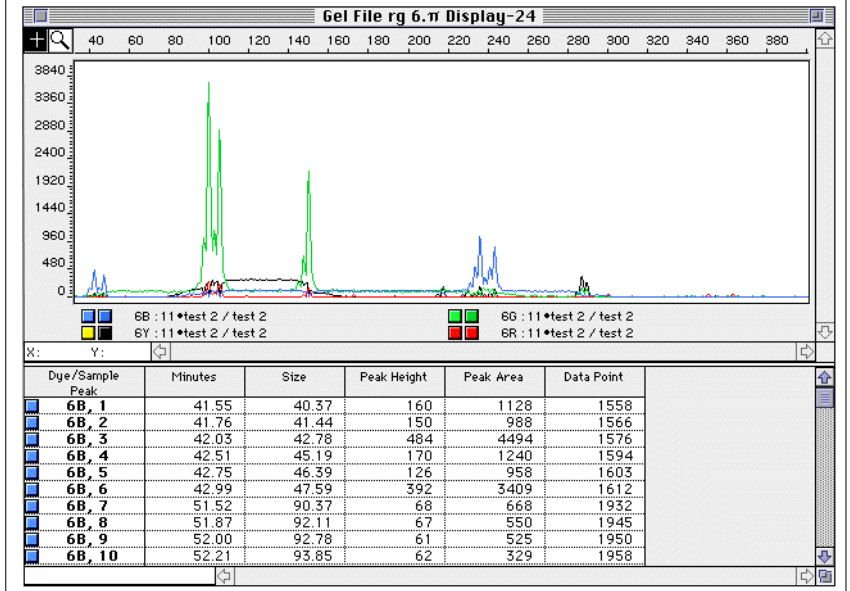
For this cause...	Take this action...
designated the wrong files.	reassign the matrix files. See "Assigning the Matrix File to Sample Files" on page 6-35.
signal is too weak to make a matrix.	rerun the matrix standards. "Loading and Running Matrix Standards on the ABI 373 with XL Upgrade" on page 6-24.

Two Causes of Bad Matrix Files The following table lists two common causes of bad matrix files:

Problem	Cause	What to do
Artifact peaks of different colors under the true peaks. See the figure below.	Loading too much dye when running matrix standards resulting in dye bleed-through.	Complete another run and recreate the matrix.



Problem	Cause	What to do
Noisy baseline.	If the matrix subtracts too much of a particular color from the sample, then the baseline may become too elevated, resulting in false peaks. See the figure below.	Complete another matrix run and make sure that you do not have any off-scale data.



Evaluating Analysis Results

7

Introduction

In This Chapter Topics in this chapter include the following:

Topic	See page
Process of Evaluating Analysis Results	7-2
Ways to Display Analysis Results	7-3
The Results Display Window	7-4
Using the Results Control Window	7-7
Changing How the Results are Displayed and Printed	7-13
The Sample Results View	7-16
Updating the Results	7-16
Saving and Renaming the Results Control Format	7-17
About Electropherogram and Tabular Data Displays	7-20
Displaying Electropherogram and Tabular Data	7-22
Displaying Electropherogram Data	7-25
Working with Electropherogram Data	7-27
How to Define Custom Colors	7-40
How to Change the Dye Scale	7-44
Process of Verifying Results	7-46
How to Verify Size Calculations	7-48
Using the Analysis Log	7-51
How to Verify Peak Detection	7-53

Process of Evaluating Analysis Results

Steps to Evaluating Analysis Results

The following table lists the steps to evaluating the analysis results:

Step	Action	For information, see...
1	Displaying Analysis Results.	"Ways to Display Analysis Results" on page 7-3.
2	Using Electropherogram and Tabular Data Displays.	"About Electropherogram and Tabular Data Displays" on page 7-20.
3	Viewing Electropherograms.	"Displaying Electropherogram Data" on page 7-25.
4	Verifying analysis results.	"Process of Verifying Results" on page 7-46.
5	Displaying Other Sample File Data.	<ul style="list-style-type: none"> ◆ "Sample Info View" on page 3-8. ◆ "Size Curve View" on page 3-13. ◆ "Raw Data View" on page 3-15. ◆ "EPT Data View" on page 3-17.
6	Using the Analysis Log.	"Using the Analysis Log" on page 7-51.
7	Remembering and Renaming the Results Display.	"Saving and Renaming the Results Control Format" on page 7-17.

Ways to Display Analysis Results

Two Ways to Display Analysis Results

The following table lists two ways to display analysis results:

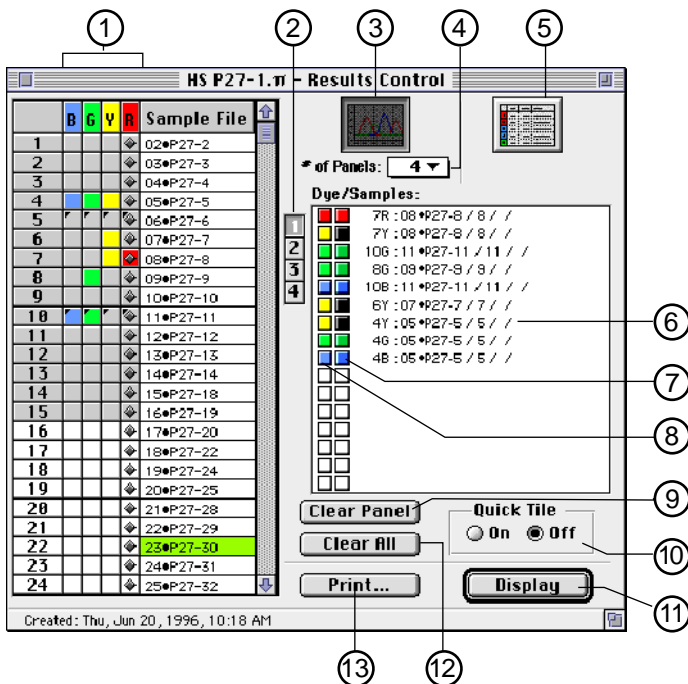
You can use...	Description	See page
The Results Display Window	This window is created from the Results Control window of a project. Use to group and view multiple Sample files as electropherogram and tabular data.	7-4
The Sample Results View	This view displays one Sample file at a time, just as the Results Display window, but is more convenient for viewing analysis results from a single Sample file. It also allows quick access to supporting information views.	7-16

The Results Display Window

Introduction The Results Display window is created from the Results Control window of a project and allows you to group and view multiple Sample files as electropherogram or tabular data. You can use the window to show up to eight panels, with multiple dye/samples per panel.

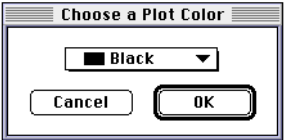
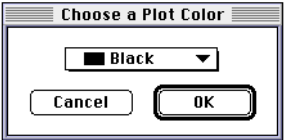
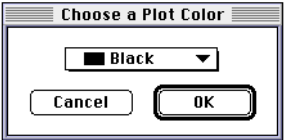
Displaying the Window Choose Results Control (⌘ 2) from the Windows menu.

The following is an example of the Results Control window.



Results Control Window Description

Results Control window descriptions:

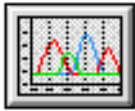

No.	Description						
1	Select dye/sample information by clicking one of the dye color fields.						
2	Each button corresponds to a panel in the display window.						
3	Click to display electropherograms for the selected samples.						
4	Choose the number of electropherogram panels available for display from the pop-up menu.						
5	Click to display tabular data for the selected samples.						
6	Identifies the sample by row number and dye code. Sample information is displayed as specified in the project options.						
7	Plot color indicator. <table border="1" data-bbox="508 613 1233 1193"> <thead> <tr> <th>If you...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>double click the plot color indicator</td> <td>the Choose a Plot Color dialog box appears.  </td> </tr> <tr> <td>⌘-double-click the dye color indicator</td> <td>the plot color indicator returns to the default color. For more information, see "Setting Dye Indicator Preferences" on page 5-15.</td> </tr> </tbody> </table>	If you...	Then...	double click the plot color indicator	the Choose a Plot Color dialog box appears. 	⌘-double-click the dye color indicator	the plot color indicator returns to the default color. For more information, see "Setting Dye Indicator Preferences" on page 5-15.
If you...	Then...						
double click the plot color indicator	the Choose a Plot Color dialog box appears. 						
⌘-double-click the dye color indicator	the plot color indicator returns to the default color. For more information, see "Setting Dye Indicator Preferences" on page 5-15.						
	Note If you change the plot color, a vertical line appears beside the indicator, showing that it has been modified.						

Results Control window descriptions:

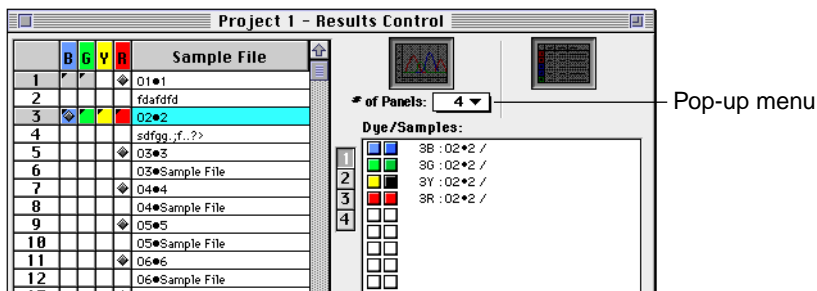
No.	Description					
8	Dye color indicator.					
	<table border="1"> <thead> <tr> <th data-bbox="452 219 821 256">If you...</th> <th data-bbox="821 219 1185 256">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="452 256 821 651">double click the dye color indicator</td> <td data-bbox="821 256 1185 651"> the Choose a Dye Scale dialog box appears. <div data-bbox="844 354 1126 493" style="text-align: center;"> </div> </td> </tr> <tr> <td data-bbox="452 651 821 821">⌘-double-click the dye color indicator</td> <td data-bbox="821 651 1185 821"> the dye color indicator returns to the default scale. <p>For more information, see “Changing the Dye Scale Preferences” on page 7-45.</p> </td> </tr> </tbody> </table>	If you...	Then...	double click the dye color indicator	the Choose a Dye Scale dialog box appears. <div data-bbox="844 354 1126 493" style="text-align: center;"> </div>	⌘-double-click the dye color indicator
If you...	Then...					
double click the dye color indicator	the Choose a Dye Scale dialog box appears. <div data-bbox="844 354 1126 493" style="text-align: center;"> </div>					
⌘-double-click the dye color indicator	the dye color indicator returns to the default scale. <p>For more information, see “Changing the Dye Scale Preferences” on page 7-45.</p>					
<p>Note If you change the scale, a vertical line appears beside the indicator, showing that it has been modified.</p>						
9	Clear panel button. See “Removing Samples” on page 7-12.					
10	Quick Tile radio buttons. See “Creating Tiled Electropherogram Displays” on page 7-10.					
11	Display button. See “Displaying the Results” on page 7-12.					
12	Clear All button. See “Removing Samples” on page 7-12.					
13	Print button. See “Printing the Results” on page 7-12.					

Using the Results Control Window

Selecting Display Format Choose the results display format by clicking either or both the electropherogram and tabular icons. You must select one of the icons to display or print data.

Click this icon...	To display...	And the...
	electropherogram data	panel information below the icon is enabled.
	tabular data	panel information below the icons is dimmed. Tabular data appears in a single table, so electropherogram panel configuration is not relevant.

Selecting Electropherogram Panels In electropherograms, the data appears in panels. You can overlay up to sixteen samples within up to eight panels. Select the number of available panels (up to eight) from the pop-up menu labeled # of Panels.



The screenshot shows the 'Project 1 - Results Control' window. On the left is a table with columns for dye (B, G, Y, R) and 'Sample File'. The table contains 12 rows of sample data. On the right is a control panel with a 'Pop-up menu' for '# of Panels' set to 4. Below this is a 'Dye/Samples' section with four color-coded buttons (B, G, Y, R) and their corresponding sample identifiers (3B, 3G, 3Y, 3R).

Note Set the number of panels when using the Quick Tile option. Using this option causes the GeneScan Analysis Software to automatically change panels as you select samples for display (see "Creating Tiled Electropherogram Displays" on page 7-10).

Selecting Samples to Display

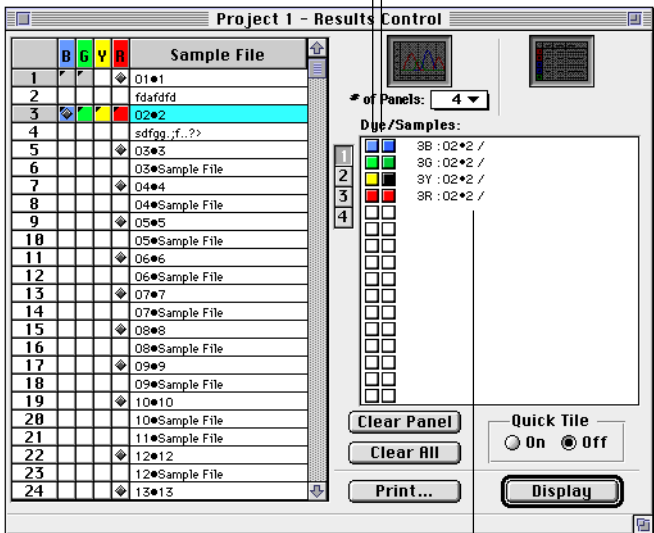
Use the dye color field on the left side of the Results Control window to select the dye for display.

You can select any or all dye colors, including the standard, for each Sample file.

To select samples for display:

Step	Action								
1	Click the Electropherogram icon. The panel below the icon is enabled.								
2	Click a panel number in the list to the left of the Dye/Sample Display list. If fewer than eight panels are available and you want to use more, choose a larger number (up to eight) from the # of Panels pop-up menu.								
3	Click dye color fields to select or deselect the corresponding samples. Take the following action: <table border="1" data-bbox="463 787 1189 982"><thead><tr><th>To...</th><th>Click the...</th></tr></thead><tbody><tr><td>select the entire column</td><td>header.</td></tr><tr><td>deselect the entire column</td><td>header, again.</td></tr><tr><td>select four colors for a Sample file</td><td>index number to the left of the color columns.</td></tr></tbody></table>	To...	Click the...	select the entire column	header.	deselect the entire column	header, again.	select four colors for a Sample file	index number to the left of the color columns.
To...	Click the...								
select the entire column	header.								
deselect the entire column	header, again.								
select four colors for a Sample file	index number to the left of the color columns.								
4	If you specified Electropherogram display, the samples corresponding to the selected dyes appear in the Dye/Sample Display list to the right of the Sample File list, as shown in the figure below.								

To select samples for display: *(continued)*

Step	Action
	<div style="display: flex; justify-content: space-around; margin-bottom: 10px;"> Dye color of sample Plot color of electropherogram </div>  <p style="text-align: right; margin-right: 50px;">Row number and dye code followed by dye/sample information</p>
5	<p>When you change to a new panel, the dye colors of the samples you selected in other panels appear dark gray to indicate that they have been selected.</p> <p>You can select them again in the current panel.</p>

Creating Tiled Electropherogram Displays

Procedure

To create tiled electropherogram displays:

Step	Action						
1	Choose the number of panels you want to display from # of Panels pop-up menu.						
2	Click the On radio button under Quick Tile.						
3	<p>Select samples by clicking color fields.</p> <p>For information on:</p> <table border="1"> <thead> <tr> <th>Topic</th> <th>See page</th> </tr> </thead> <tbody> <tr> <td>Selecting Samples to Display</td> <td>7-8</td> </tr> <tr> <td>Setting the tiled electropherogram preferences</td> <td>7-13</td> </tr> </tbody> </table> <p>Each time a sample is selected, the program automatically changes to the next panel, so each selection is placed after the one containing your previous selection.</p> <p>After you select a sample for the last panel, the panel increments to the first panel again.</p>	Topic	See page	Selecting Samples to Display	7-8	Setting the tiled electropherogram preferences	7-13
Topic	See page						
Selecting Samples to Display	7-8						
Setting the tiled electropherogram preferences	7-13						

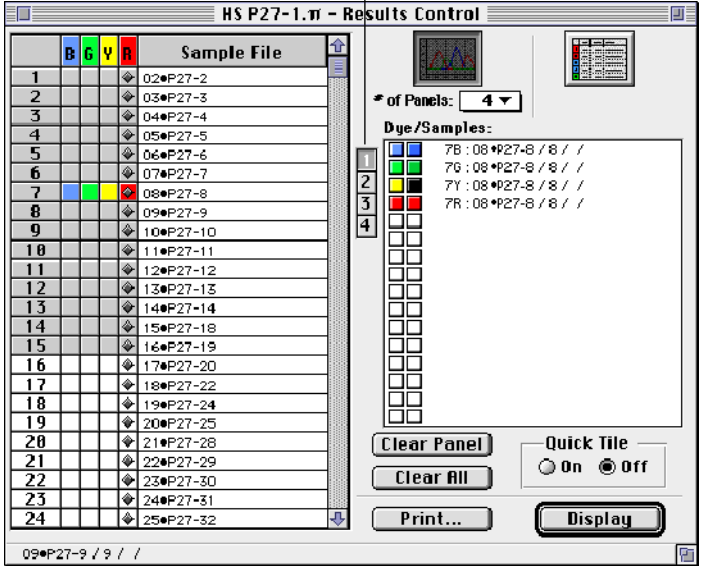
Example

The following table describes two examples of how to use the Quick Tile feature:

If you...	Then...
have four samples and choose four panels for display	<p>click the column heading for the blue dye to select all four blue dye-labeled samples.</p> <p>The blue dye for each Sample file appears in a separate panel.</p>
click the row index number for the first Sample file to select all dye colors for one sample file	each dye for that Sample file appears in a separate panel.

Deselecting Samples

To deselect the samples that you have selected for display:

Step	Action
1	<p>If you specified Electropherogram display, click the panel number to the left of the Dye/Samples display list to display the information for the panel containing the samples you want to deselect.</p> <p style="text-align: center;">Dye color fields Panel numbers</p>  <p>The screenshot shows the 'HS P27-1.n - Results Control' window. On the left, there is a table with columns for dye colors (B, G, Y, R) and 'Sample File'. The table lists samples from 02 to 25. Sample 08 is highlighted with a yellow background. To the right of the table is a 'Dye/Samples' display area with a grid of colored squares and a list of sample identifiers. Below the grid are buttons for 'Clear Panel', 'Clear All', 'Print...', and 'Display'. A 'Quick Tile' section has 'On' and 'Off' radio buttons.</p>
2	<p>On the left side of the Results Control window, deselect the dye color fields corresponding to the samples you want to remove.</p> <p>The dye/sample identifiers are removed from the Dye/Sample Display list as you deselect samples.</p>

Removing Samples

To remove samples:

Step	Action						
1	Select the panel number on the buttons to the left of the Dye/Sample Display list.						
2	You can take the following action: <table border="1"><thead><tr><th>To remove...</th><th>Click...</th></tr></thead><tbody><tr><td>all the samples you selected for display in a panel</td><td>Clear Panel.</td></tr><tr><td>the samples you have selected for display in all panels</td><td>Clear All.</td></tr></tbody></table>	To remove...	Click...	all the samples you selected for display in a panel	Clear Panel.	the samples you have selected for display in all panels	Clear All.
To remove...	Click...						
all the samples you selected for display in a panel	Clear Panel.						
the samples you have selected for display in all panels	Clear All.						

Displaying the Results

To display the results on the screen, you can either:

- ◆ Click the Display button.
- ◆ Press the Return or Enter key.

Printing the Results

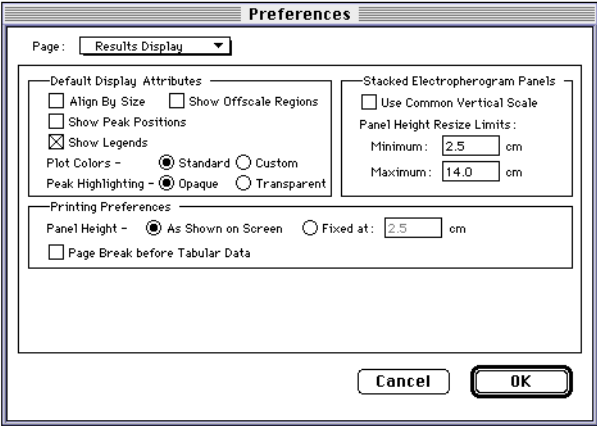
To print the results:

Step	Action
1	You can either: <ul style="list-style-type: none">◆ Click the Print button, or◆ Choose Print (⌘ P) from the File menu.
2	Click Print in the dialog box that appears.

Changing How the Results are Displayed and Printed

Procedure You can set certain display preferences that remain in effect each time you display or print results data.

To change how results are displayed and printed:

Step	Action
1	<p>Choose Preferences from the Settings menu and choose Results Display from the submenu.</p> <p>The following dialog box appears.</p>  <p>There are three Results Display preferences:</p> <ul style="list-style-type: none">◆ Default Display Attributes.◆ Stacked Electropherogram Panels.◆ Printing Preferences.

To change how results are displayed and printed: *(continued)*

Step	Action	
<p>2</p>	<p>Set the Default Display Attributes to control the display attributes of new results displays, as follows:</p>	
	<p>You can set...</p>	<p>For more information...</p>
	<p>Show Peak Positions</p>	<p>“Displaying Peak Positions” on page 7-29.</p>
	<p>Use opaque or transparent peak highlighting</p>	<p>“Highlighting Peaks” on page 7-29.</p>
	<p>Show Legends</p>	<p>“Using Legends to Change the Display” on page 7-30.</p>
	<p>Show Offscale Data</p>	<p>“Showing Off-Scale Data” on page 7-32.</p>
	<p>Align By Size</p>	<p>“Showing Data by Fragment Size” on page 7-35.</p>
	<p>Use standard or custom plot colors</p>	<p>“Defining Custom Color for All Electropherogram” on page 7-41.</p>
<p>3</p>	<p>Configure Stacked Electropherogram Panels, as follows:</p>	
	<p>Choose...</p>	<p>To...</p>
	<p>Use Common Vertical Scale checkbox</p>	<p>set all panels in a display so they have the same vertical scale.</p> <p>The common scale is based on the electropherogram with the largest vertical scale.</p>
<p>Panel Height Resize Limits</p>	<p>set minimum and maximum values for electropherogram panel height in the results display.</p> <p>This allows you to limit how much the electropherogram panels stretch or shrink to fit the size of the window.</p>	

To change how results are displayed and printed: *(continued)*

Step	Action		
4	Set the Printing Preferences, as follows:		
	<table border="1"><thead><tr><th data-bbox="512 250 874 289">If...</th><th data-bbox="874 250 1245 289">Then...</th></tr></thead></table>	If...	Then...
	If...	Then...	
	the height of the panels that appears on screen is acceptable	choose the As Shown on Screen radio button.	
you want to print the electropherogram at a specified height	click the Fixed at radio button and enter a value in the field.		
you want to force a page break after the electropherograms have been printed	select the Page Break before Tabular data checkbox.		
5	Click OK		

The Sample Results View

About the View The Sample Results view is displayed within a Sample file window. You can access the window through a Sample file or through a project's Analysis Control or Results Control window. If you are opening a Sample file as a stand-alone file, the Sample Results view is the default display within the Sample file window.

For More Information See "Sample Results View" on page 3-6.

Updating the Results

Re-Analyzing the Data The Results Control and the Sample Results windows are dynamic. The following table lists how the windows are updated if you re-analyze the data.

If you re-analyze...	Then the...
your data with either window active	software updates the window.

Saving and Renaming the Results Control Format

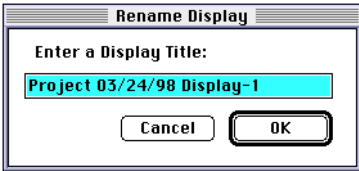
Introduction You can use the Results Control window to view multiple Sample files in electropherogram and tabular format. The GeneScan Analysis Software allows you to save formats for future use. You can then redisplay or print these formats without having to redefine them again.

Important Considerations The following are important considerations for saving a Results Control format:

- ◆ You must save the project for the display to be available when you open the project again.
- ◆ Remembering a display preserves the combination of windows/panels /data and customized color settings. It does not preserve any zooming you have performed.

Saving the Display Format

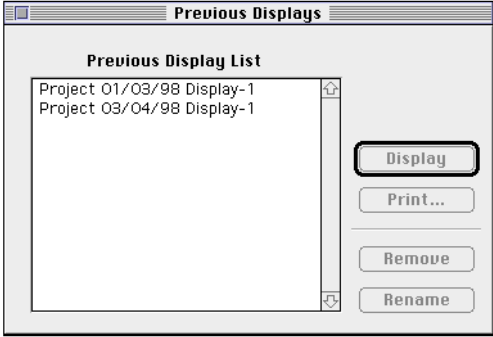
To save a display format for future viewing:

Step	Action
1	With the Results Control window set for the display, either: <ul style="list-style-type: none">◆ Click the Display button, or◆ Press the Return key.
2	With the display on the screen, choose Remember Display from the Project menu. The Remember Display dialog box appears. 
3	Enter a name for the display and click OK.

Working with a Previously Saved Displays

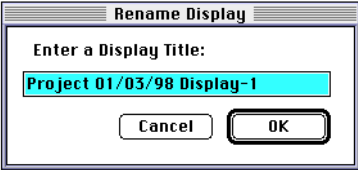
Use the Previous Displays dialog box to display, print, remove or rename a saved display.

To work with a previously saved displays:

Step	Action										
1	<p>Choose Previous Displays from the Project menu.</p> <p>The following is an example of the Previous Displays dialog box.</p> 										
2	<p>Select a display or multiple displays and take the following action:</p> <table border="1" data-bbox="467 824 1185 1380"> <thead> <tr> <th data-bbox="467 824 826 894">To...</th> <th data-bbox="826 824 1185 894">Click an item in the list and click...</th> </tr> </thead> <tbody> <tr> <td data-bbox="467 894 826 935">display the saved formats</td> <td data-bbox="826 894 1185 935">Display.</td> </tr> <tr> <td data-bbox="467 935 826 1049">print the saved formats</td> <td data-bbox="826 935 1185 1049">Print. The standard print dialog box appears.</td> </tr> <tr> <td data-bbox="467 1049 826 1130">remove the saved formats</td> <td data-bbox="826 1049 1185 1130">Remove. An alert appears.</td> </tr> <tr> <td data-bbox="467 1130 826 1380">rename</td> <td data-bbox="826 1130 1185 1380">Rename. Note You can only rename one display at a time. The Rename dialog box appears (See “Renaming the Current Results Display” on page 7-19).</td> </tr> </tbody> </table>	To...	Click an item in the list and click...	display the saved formats	Display.	print the saved formats	Print. The standard print dialog box appears.	remove the saved formats	Remove. An alert appears.	rename	Rename. Note You can only rename one display at a time. The Rename dialog box appears (See “Renaming the Current Results Display” on page 7-19).
To...	Click an item in the list and click...										
display the saved formats	Display.										
print the saved formats	Print. The standard print dialog box appears.										
remove the saved formats	Remove. An alert appears.										
rename	Rename. Note You can only rename one display at a time. The Rename dialog box appears (See “Renaming the Current Results Display” on page 7-19).										

Renaming the Current Results Display

To rename the Results display that is currently on the screen and to save the display under a different name.

Step	Action
1	Ensure the display is the active window.
2	Choose Rename Display from the Project menu. The following is an example of the Rename Display dialog box. 
3	Enter a new name for the display and click OK.

About Electropherogram and Tabular Data Displays

Introduction After analyzing the data, you can display the results for each sample in electropherograms and tabular data. You can also customize the electropherograms and tabular data display.

Note Altering the appearance of the electropherograms and the tabular data displays does not change the analyzed data contained in the Sample file on which they are based.

How the Window is Divided When electropherograms and tabular data are displayed together, the window is divided into upper and lower windows.

Window	Contains
Upper window	electropherograms
Lower window	tabular data

You can customize the window's appearance by adjusting the relative size of each window, see "Adjusting Window Size" on page 7-24.

What Tabular Data Contains If you analyze samples with an internal size standard and perform size calling, the tabular data contains the estimated sizes (in base pairs) of all detected fragments. Use this information for detailed data analysis and further calculations. The peaks matched to the defined size standard are identified by dots next to the Dye/Sample Peak field.

Sample peaks that are larger (in base pairs) than the largest defined peak in the selected standard are not sized. The corresponding size fields are blank.

Note Tabular data displays only peaks that are detected based on the Dye Amplitude Thresholds and Minimum Peak Half Width setting of the analysis parameters.

How Electropherogram Panels are Sized In the Results Display window, the GeneScan Analysis Software sizes all electropherogram panels to fit within the electropherogram portion of the window by using the largest size that fits them into the visible area. You can scroll to see the portion of the display that is not visible.

**Peaks Visible in
Electropherogram
Not Tabular Data**

Peaks may be visible in the electropherogram and not listed in the tabular data because:

Reason	For more information, see
The software detects the peaks based on the Dye Amplitude Thresholds and Min Peak Half Width.	“Peak Detection Parameter Options” on page 5-20.
Electropherograms display the peaks that fall within the range specified by the Size Call Range parameters that are defined in the Analysis Parameters dialog box.	“Size Call Range Parameter Options” on page 5-21.

**For More
Information**

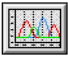

For more information, see the following topics:

Topic	See
Displaying Electropherogram and Tabular Data	7-22
Displaying Electropherogram Data	7-25
Working with Electropherogram Data	7-27

Displaying Electropherogram and Tabular Data

Displaying Electropherogram and Tabular Data

To display electropherogram and tabular data:

Step	Action
1	Choose Results Control (☼ 2) from the Windows menu to open the Results Control window.
2	Click the electropherogram icon () and the tabular icon ()
3	If applicable, select the number of electropherogram panels.
4	Select the dye/samples for display.
5	Click Display. The electropherogram and tabular data are displayed in the Results Display window. See also “Working with Electropherogram Data” on page 7-27.

Tabular Data and Electropherogram Example

The following is an example of tabular data with a corresponding electropherogram:

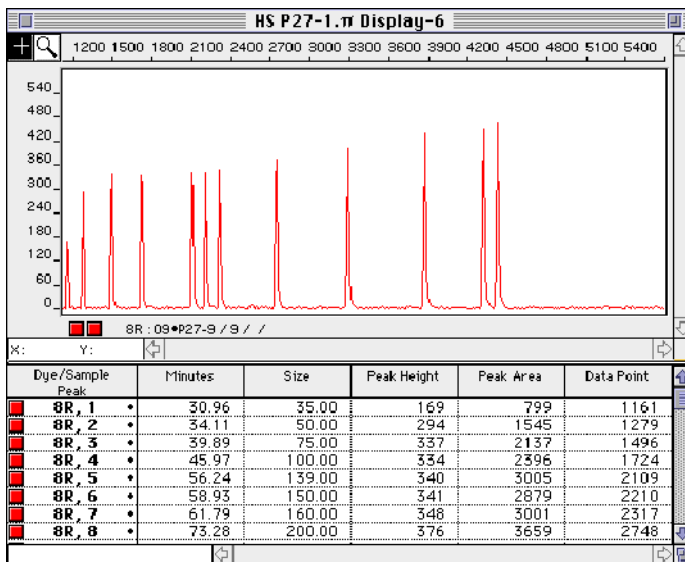


Table Describing Columns

The following table describes the columns in the above figure:

This column	Identifies...
Dye/Sample Peak	<ul style="list-style-type: none">◆ Sample index number.◆ Dye color.◆ Peak number.
Minutes	The time, in minutes, from the start of the run to the time the fragment was detected.
Size	The differences in fragment mobility. This value is calculated automatically only if you: <ul style="list-style-type: none">◆ Run the size standard in the same lane or injection as the sample, and◆ Perform size calling.
Peak Height	Signal size.
Peak Area	Area of the detected peak.
Data point	Data point of the fragment at its maximum peak height.

Highlighting Information

To highlight information for one peak in the electropherogram and tabular data:

Click...	Then...
the peak in the electropherogram	the peak fills with color and the corresponding row in the tabular data window is highlighted.
the Dye/Sample Peak number in the tabular data window	highlights the corresponding peak in the electropherogram.


How to Change the Highlight Color

Choose Peak Highlighting from the View menu and either Opaque or Transparent from the submenu.

For more information, see “Highlighting Peaks” on page 7-29.

Adjusting Window Size

To adjust the relative size of the electropherogram and tabular windows:

Step	Action
1	Move the cursor to the window divider (the double line between the two windows).
2	When the cursor changes to a bidirectional arrow (), click the window divider line and drag it up or down.

Hiding Selected Rows of Data

To hide selected rows of data:

Step	Action								
1	Take the following action: <table border="1"><thead><tr><th>If you want to...</th><th>Then...</th></tr></thead><tbody><tr><td>select a row</td><td>either:<ul style="list-style-type: none">◆ Click the first field in the row, or◆ Click the corresponding peak in the electropherogram.</td></tr><tr><td>select several consecutive rows</td><td>shift-click the first and last row you want to select.</td></tr><tr><td>select several rows that are not next to each other</td><td>⌘-click the rows.</td></tr></tbody></table>	If you want to...	Then...	select a row	either: <ul style="list-style-type: none">◆ Click the first field in the row, or◆ Click the corresponding peak in the electropherogram.	select several consecutive rows	shift-click the first and last row you want to select.	select several rows that are not next to each other	⌘-click the rows.
If you want to...	Then...								
select a row	either: <ul style="list-style-type: none">◆ Click the first field in the row, or◆ Click the corresponding peak in the electropherogram.								
select several consecutive rows	shift-click the first and last row you want to select.								
select several rows that are not next to each other	⌘-click the rows.								
2	Choose Hide Selected Rows (⌘ H) from the View menu.								

Limited the Rows Displayed

To limit the display to the selected rows of data:

Step	Action
1	Select the rows you want to display.
2	Choose Show ONLY selected Rows (⌘ G) from the View menu. Note Choose Show All Rows (⌘ G) from the View menu to display all of the tabular data after limiting the display.

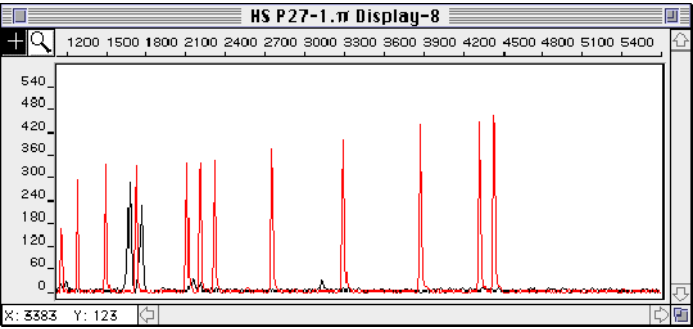
Displaying Electropherogram Data

Definition Each electropherogram provides a profile of the selected dye samples it represents. The y-axis represents the relative fluorescence of the detected fragments as they occurred over time. The x-axis represents time and can be displayed by data points or base pairs.

Base Pairs Versus Data Points You can choose to have the horizontal tick marks on the x-axis represent size in base pairs instead of data points only if you ran an internal size standard with the sample (see “Showing Data by Fragment Size” on page 7-35).

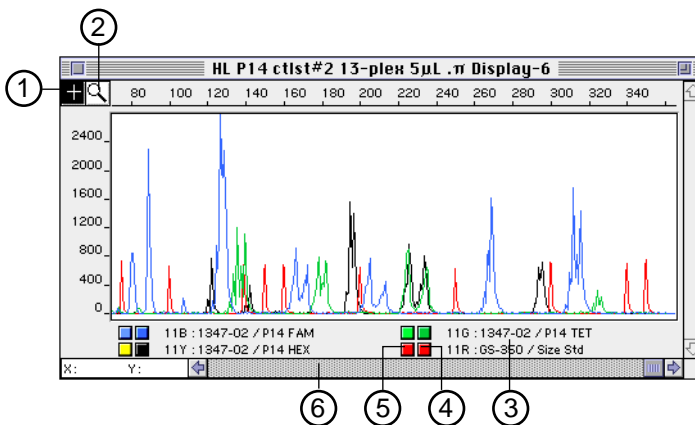
Procedure To display electropherogram data:

Step	Action
1	Choose Results Control (☞ 2) from the Windows menu to open the Results Control window.
2	Click the electropherogram icon.
3	If applicable, select the number of electropherogram panels.
4	Select the dye/samples for display.
5	Click Display. The electropherogram is displayed in the Results Display window.



Electropherogram Example

The following is an example of an electropherogram:



Descriptions of the Electropherogram

The following table lists the descriptions for the above figure:

No.	Description	See...
1	Cross hairs	“Displaying X-and Y-Axis Positions” on page 7-27.
2	Magnifying glass Use to zoom in a specific area or hold down while pressing the Option key to zoom out in successively smaller scale.	“Zooming In and Out” on page 7-31.
3	Legend Text from the Sample file that appear beneath electropherogram panels in the Results Display window.	“Using Legends to Change the Display” on page 7-30.
4	Dye color indicator	“Dye color indicator.” on page 7-6.
5	Plot color indicator	“Plot color indicator.” on page 7-5.
6	Scroll bar Use the scroll bar to scroll horizontally.	“Scrolling the Display” on page 7-30.

Working with Electropherogram Data

In This Section This section describes how to perform the following tasks:

Task	See
Displaying X-and Y-Axis Positions	7-27
Moving the Electropherogram	7-28
Changing the Dye Color	7-28
Displaying Peak Positions	7-29
Highlighting Peaks	7-29
Using Legends to Change the Display	7-30
Scrolling the Display	7-30
Zooming In and Out	7-31
Showing Off-Scale Data	7-32
Showing Data by Fragment Size	7-35
Changing the Horizontal Scale	7-37
Changing the Vertical Scale	7-38
Assigning Standard or Custom Colors	7-39

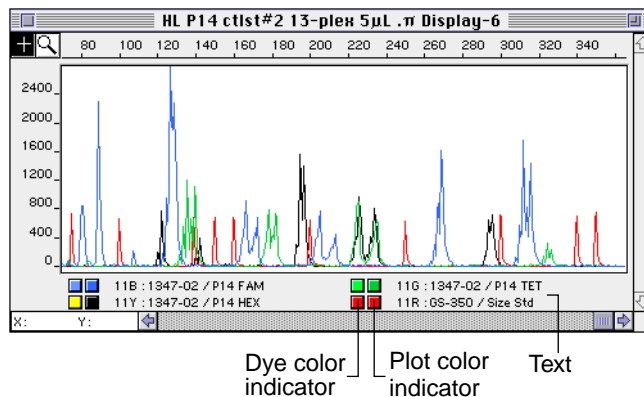
Displaying X-and Y-Axis Positions Click the cross hairs and select an area in the Electropherogram to display the x-and y-axis positions.

The x-and y-axis values appear in the box in the lower left corner of the electropherogram and, if tabular data is also displayed, the row in the table is highlighted.

Moving the Electropherogram

Move the associated electropherogram to the front by clicking in the legend either the:

- ◆ Dye color indicator
- ◆ Plot color indicator, or
- ◆ Text



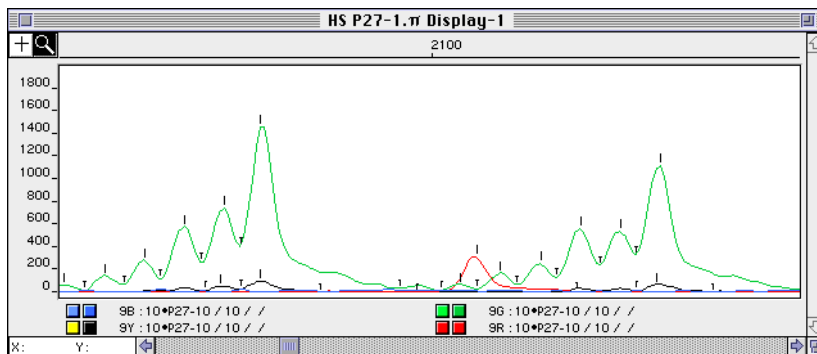
Changing the Dye Color

To change the dye color, double-click the dye color indicator. If you change the dye color or scale, a vertical line appears beside the indicator, showing that it has been modified.

Note ☞-double-click the dye color indicator to return it to the default.

Displaying Peak Positions

Choose Show Peak Positions from the View menu. Use this command to examine how the GeneScan Analysis Software defined peaks by displaying markers that identify the beginning, center, and end of each peak.



Highlighting Peaks

Use the Peak Highlighting command to highlight a selected peak with the dye/sample's plot color.

To highlight selected peaks:

Step	Action						
1	Choose Peak Highlight from the Views menus and either Opaque or Transparent from the submenu.						
	<table border="1"> <thead> <tr> <th>Use this option...</th> <th>To...</th> </tr> </thead> <tbody> <tr> <td>Opaque</td> <td>fill the peak with a solid color that can obscure peaks behind the selected peak.</td> </tr> <tr> <td>Transparent</td> <td>use a slightly diffused plot color that allows you to view overlapping peaks.</td> </tr> </tbody> </table>	Use this option...	To...	Opaque	fill the peak with a solid color that can obscure peaks behind the selected peak.	Transparent	use a slightly diffused plot color that allows you to view overlapping peaks.
	Use this option...	To...					
Opaque	fill the peak with a solid color that can obscure peaks behind the selected peak.						
Transparent	use a slightly diffused plot color that allows you to view overlapping peaks.						
2	Click a detected peak in an electropherogram. The peak is highlighted						



Using Legends to Change the Display

The following table lists how to use legends to change how electropherograms are displayed:

If you want to...	Then...
show or hide legends	choose Show Legends from the View menu.
open Sample file windows	double-click the corresponding legend text.
reorganize overlaid electropherograms	<ol style="list-style-type: none"> Display the electropherograms with legends. Click either the dye, color indicator, plot color indicator, or the text for the sample you want to move to the front.

Scrolling the Display

The following table lists ways to scroll the display:

Use the	Description								
scroll bar	<table border="1"> <thead> <tr> <th>If you want to...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>shift the electropherogram to the right or the left</td> <td>click in the gray region of the scroll bar to the right or the left of the scroll box.</td> </tr> <tr> <td>scroll across the electropherogram</td> <td>click an arrow at the end of the scroll bar.</td> </tr> <tr> <td>control the amount of scroll</td> <td>drag the scroll bar to the right or the left.</td> </tr> </tbody> </table>	If you want to...	Then...	shift the electropherogram to the right or the left	click in the gray region of the scroll bar to the right or the left of the scroll box.	scroll across the electropherogram	click an arrow at the end of the scroll bar.	control the amount of scroll	drag the scroll bar to the right or the left.
	If you want to...	Then...							
	shift the electropherogram to the right or the left	click in the gray region of the scroll bar to the right or the left of the scroll box.							
	scroll across the electropherogram	click an arrow at the end of the scroll bar.							
control the amount of scroll	drag the scroll bar to the right or the left.								
scroller symbols	<ol style="list-style-type: none"> Hold the mouse cursor over either the vertical or horizontal scale of the electropherogram. Either a vertical scroller symbol () or a horizontal scroller symbol () appears. Hold down the mouse button and move the mouse in the direction of the information you want to view. 								

Zooming In and Out

Out

About Zooming In and Out

By default the GeneScan Analysis Software scales each electropherogram horizontally to show all peaks detected during the run. While this provides a good overview of the run, some peaks may be quite compressed.

To improve visibility, you can change the horizontal scale of the electropherograms by zooming.

Zooming affects:

- ◆ Only the horizontal scale, and zooms the middle portion of the window.
- ◆ All displayed electropherogram panels.

How to Change the View Scale

To change the view scale:

If you want to...	Then...
see views with greater detail	<ul style="list-style-type: none">◆ Choose Zoom In (⌘ +) from the View menu, or◆ Click the magnifying glass in the upper-left corner of the window and select a specific area on which to zoom in by drawing a box on the electropherogram.
see a smaller scale view of the data after zooming in	<ul style="list-style-type: none">◆ Choose Zoom Out (⌘ -) from the View menu, or◆ Click the magnifying glass cursor, hold down the Option key, and click the electropherogram. The data appears in successively smaller scale views.
quickly scale the data so that the entire length fits within the window, again	Choose Zoom Out (Full Range) from the View menu.

Showing Off-Scale Data

This section contains the following information:

Topic	See page
Procedure	7-32
For ABI PRISM 377 Runs	7-33
About Flat Topped Peaks	7-33
Electropherogram Displaying Off-Scale Data	7-34
Electropherogram Displaying the Flat Topped Effect	7-35

Procedure

To show off-scale data:

Step	Action
1	<p>Choose Preferences from the Settings menu and Results Display from the submenu.</p> <p>The Results Display Preferences dialog box appears.</p>
2	<p>Choose the Show Offscale Regions checkbox to highlight with a red bar regions in the electropherogram that contain off-scale (see “Electropherogram Displaying Off-Scale Data” on page 7-34).</p> <p>Note Choose the Zoom In command from the View menu to more clearly show the areas of off-scale data.</p> <p>The Analysis Log lists the numbers of off-scale regions in the analysis range for each Sample file, if the sample was sized.</p> <p>Note You can toggle this command for individual electropherograms by choosing Hide/Show Offscale regions from the View menu.</p>

For ABI PRISM 377 Runs

The Show Offscale Regions feature must be used in conjunction with:

For ABI PRISM 377 runs, use	Refer to...
Pre-averaging offscale detection in the Gel Preferences menu	"Pre-Averaging Offscale Detection checkbox" on page 2-9
Suppress Left/Right Averaging checkbox in the General Setting Preferences dialog box	ABI PRISM 377 Data Collection software. Refer to Chapter 4 in the ABI PRISM 377 Instrument User's Manual.

About Flat Topped Peaks

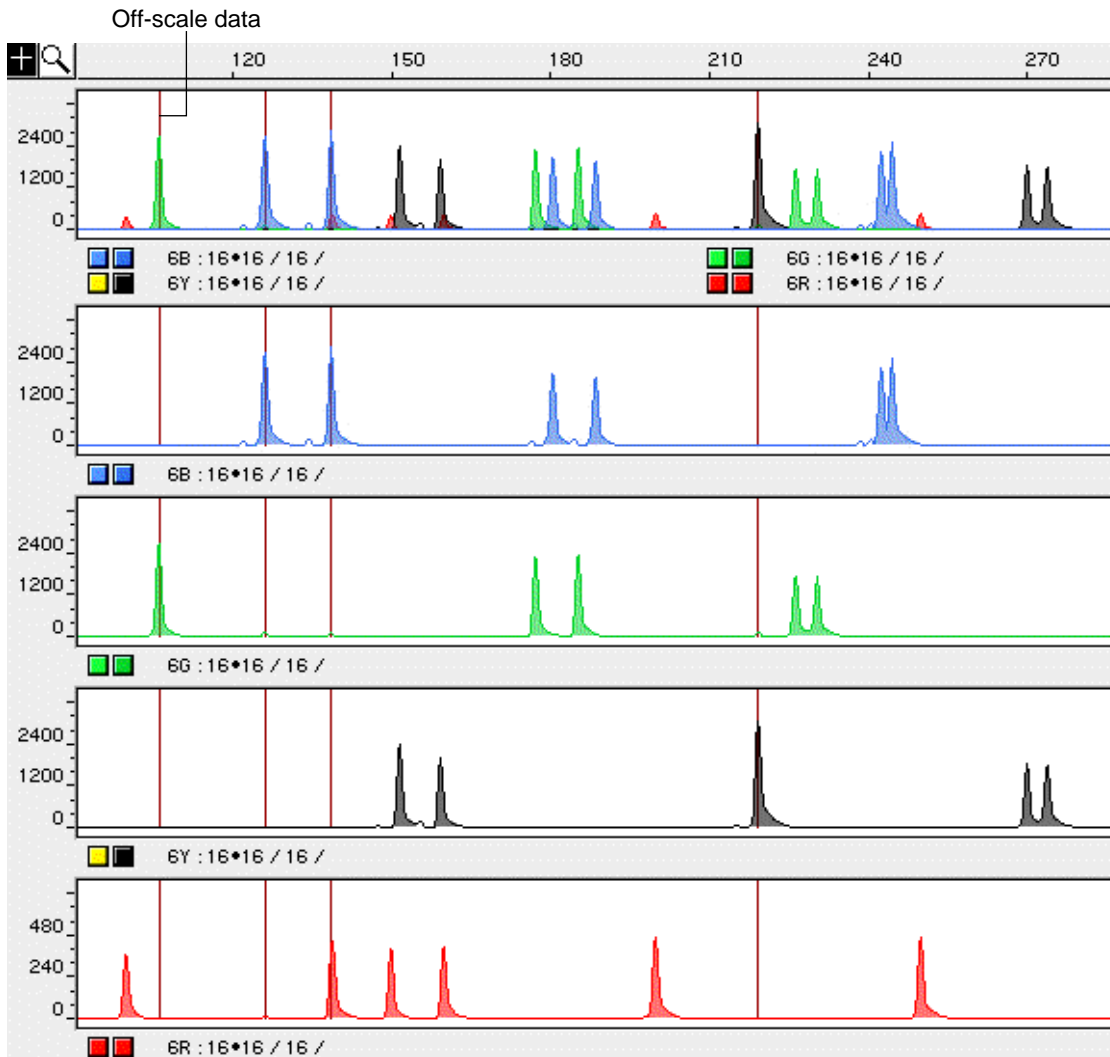
An additional feature is that peaks that contain off-scale data points are drawn in the electropherograms as "flat topped," that is, the top section of the peak is flat rather than pointed (see "Electropherogram Displaying the Flat Topped Effect" on page 7-35).

This feature may be seen when the data is analyzed with no or light smoothing; the flat topped peaks may not be apparent with heavy smoothing.

Choose Analysis Parameters from the Settings Menu (see "Smooth Options" on page 5-20).

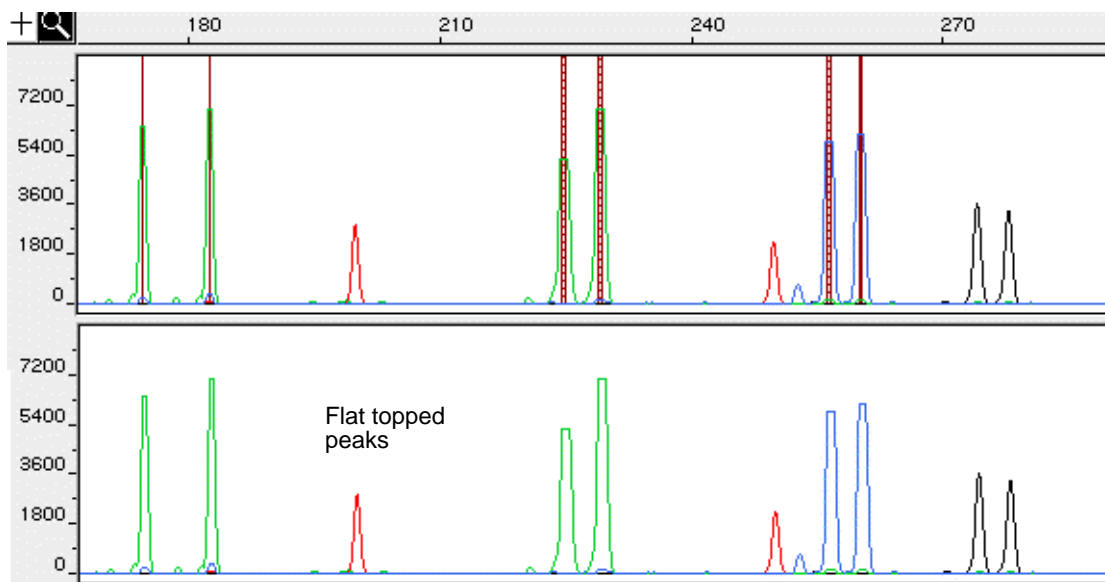
Electropherogram Displaying Off-Scale Data

The following is an example of an electropherogram displaying off-scale data.



Electropherogram Displaying the Flat Topped Effect

The off-scale peaks in the electropherogram in the expanded view below illustrates the flat topped effect.



Showing Data by Fragment Size

Introduction

If you analyze your samples with an internal size standard, you can use the Align by Size command to show the horizontal scale of the electropherograms by fragment size instead of by data point.

Note You can only display data by size if you analyzed your samples using an internal size standard.

Adjusts for Run Variations

The Align by Size command adjusts for run to run and lane to lane variations and thereby eliminates any apparent differences, and any possible confusion, caused by run discrepancies. For example, if you run two identical samples in different runs, aligning peaks by size value eliminates any apparent differences (and any possible confusion) caused by run discrepancies. With ABI 373 and ABI PRISM 377 data, the command lets you align two different samples run on the same gel using the internal size standard as a reference.

Note You can display overlaid samples in the same dye in different colors. See “How to Define Custom Colors” on page 7-40.

How to Switch Between Size and Data Point Display

To switch between size and data point display:

If you want to...	Then...
show data by size	choose Align by Size (≡ T) from the View menu. When the data is aligned by size, the menu changes to Align by Data point. Choose the command again, to show the data aligned by data point.
set the default peak alignment	choose Preferences from the Settings menu and choose Results Display from the submenu. You can use the Results Display Preferences dialog box to set certain preferences that remain in effect each time you display or print results data. For more information, see “Changing How the Results are Displayed and Printed” on page 7-13.

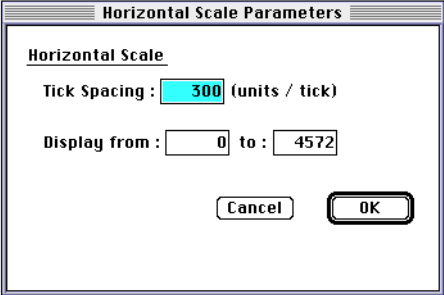
Changing the Horizontal Scale

You can change the vertical scale:

- ◆ All electropherograms.
- ◆ Individual electropherograms.

How to Change the Horizontal Scale for All Electropherograms

To change the scale of the horizontal axis for all electropherograms:

1	Display the electropherogram panels you want to change.
2	<p>Choose Horizontal Scale from the View menu.</p> <p>The Horizontal Scale Parameters dialog box appears.</p> <p>Note You can also move the cursor over the horizontal axis of a displayed electropherogram and double-click.</p> 
3	Enter the increments represented by the tick marks for the horizontal axis in the Tick Spacing box.
4	Enter a range in the entry fields labeled Display from and to.
5	Click OK.

How to Change the Horizontal Scale for Individual Electropherograms

To change the horizontal scale for individual electropherograms:

Step	Action
1	Display the electropherograms.
2	Move the cursor over the horizontal axis of the panel that you want to change and double-click. The Horizontal Scale Parameters dialog box appears.
3	Enter the increments represented by the tick marks for the horizontal axis in the Tick Spacing box.
4	Enter a range in the entry fields labeled Display from and to.
5	Click OK. The horizontal scale changes

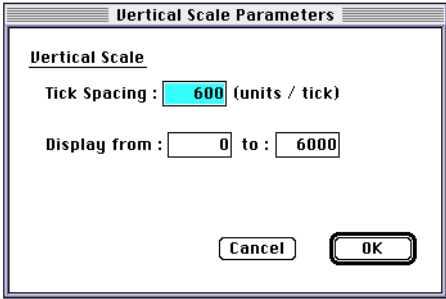
Changing the Vertical Scale

You can change the vertical scale for

- ◆ All electropherograms.
- ◆ Individual electropherograms.

How to Change the Vertical Scale for All Electropherograms

To change the vertical scale for all electropherograms:

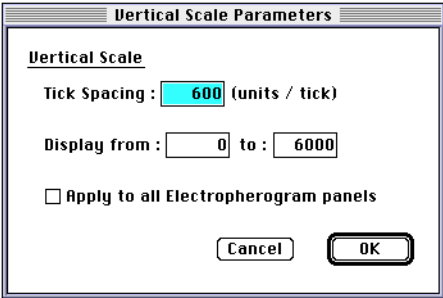
Step	Action
1	Display the electropherogram panel that you want to change.
2	Choose Vertical Scale from the View menu. The Vertical Scale dialog box appears. 
3	Enter the increments represented by the tick marks for the vertical axis in the Tick Spacing box.

To change the vertical scale for all electropherograms: *(continued)*

Step	Action
4	Enter a range in the entry fields labeled Display from and to.
5	Click OK.

How to Change the Vertical Scale for Individual Electropherograms

To change the vertical scale for individual electropherograms.

Step	Action
1	Display the electropherograms.
2	Move the cursor over the vertical axis of the panel that you want to change and double-click. The following dialog box appears.
	
3	Enter tick mark increments and a range.
4	Ensure that the Apply to all Electropherogram panels checkbox is not selected. Select the checkbox only to apply the changes to all displayed electropherogram panels.
5	Click OK. The vertical scale changes only for the electropherogram panel that you selected.

Assigning Standard or Custom Colors

To assign standard or custom colors, choose the Plot Color command from the View menu and either Standard or Custom from the submenu.

For information on defining custom colors, see “How to Define Custom Colors” on page 7-40.

How to Define Custom Colors

Introduction The GeneScan Analysis Software assigns a plot color to each dye/sample added to an electropherogram. Normally, it is the color associated with the individual dye/sample by the Dye Indicators Preferences.

Note To change the default dye colors in the Analysis Control window and the Results displays, refer to “Setting Dye Indicator Preferences” on page 5-15.

Note Custom plot colors are not available in the Sample Results view.

Why Change Colors in the Electropherogram

Change the colors in the electropherogram to:

- ◆ Differentiate between different samples labeled with the same color dye.
 - ◆ Improve contrast between different dye colors.
 - ◆ Show data in a special color for a presentation.
 - ◆ Optimize plot colors for a particular printer.
-

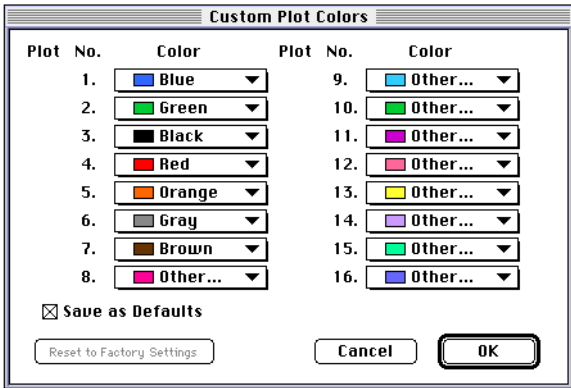
Saving the Display Format

The following table lists the options for saving the display format after customizing the display colors:

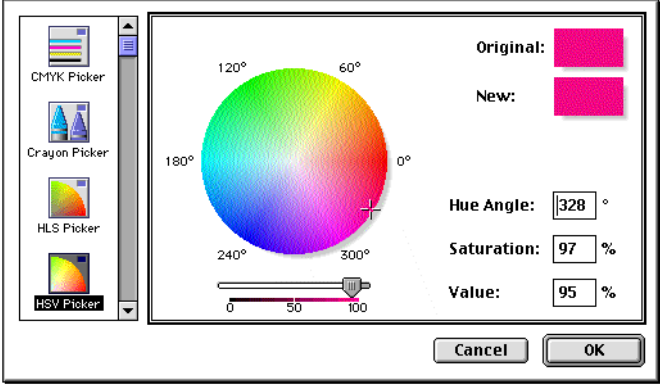
If you...	And then...
specify save the display format of a Results Display window after customizing the display	open it at a later time, the custom colors still appear. For more information, see “Saving and Renaming the Results Control Format” on page 7-17.
do not save the display format after manually customizing the colors	redisplay the same results again, the electropherograms are redrawn with default colors.

Defining Custom Color for All Electropherogram

To define custom colors for all electropherograms:

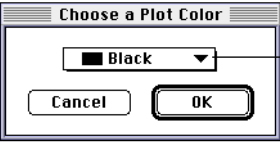
Step	Action
1	<p>Choose Project Options from the Settings menu and Choose Custom Plot Colors from the submenu.</p> <p>The Custom Plot Colors dialog box appears.</p> 
2	<p>Choose new colors from the pop-up menus beside the plot numbers.</p> <p>Note The plot numbers indicate the order of the samples in the electropherogram legend.</p>

To define custom colors for all electropherograms: *(continued)*

Step	Action
3	<p>Choose Other from the pop-up menu to specify a color that does not appear in the pop-up menu.</p> <p>The following is an example of a color picker that appears.</p> 
4	Select the checkbox labeled Save As Defaults to save the customized colors.
5	Click OK.

Defining Individual Plot Colors

To define the individual plot colors:

Step	Action
1	Display the electropherograms with legends.
2	<p>Double-click the plot color indicator next to the sample you want to change.</p> <p>The Choose a Plot Color dialog box appears.</p> 
3	<p>Choose a new color from the pop-up menu.</p> <p>If you choose Other, then a color picker appears.</p>

Step	Action
4	<p data-bbox="521 142 619 167">Click OK.</p> <p data-bbox="521 188 1224 269">The color of the electropherogram for the individual sample changes, and a vertical line appears beside the plot color indicator to signify that it is modified.</p> <p data-bbox="521 293 1224 407">Note You can change the plot color in the same way from the Results Control window. When you do so, the dye/sample is plotted with the set color each time you open the applicable Results Display window.</p> <p data-bbox="521 431 1224 480">Note Press ⌘ and double-click the plot color indicator to reset it to the original color.</p>

How to Change the Dye Scale

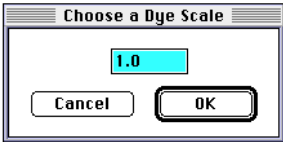
What the Dye Scale Defines The dye scale defines how dyes in an electropherogram appear relative to each other. You can compensate for peaks with different intensities by redefining the dye scale.

Note Changing the dye scale affects only the display, not the underlying data.

Increasing the Dye Scale Example The following table describes an example of increasing the dye scale.

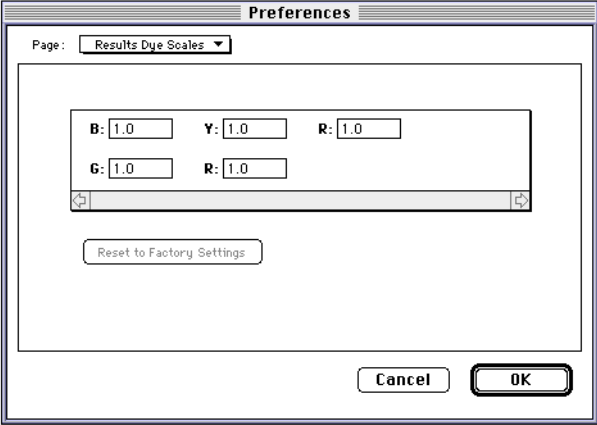
If...	Then...	Action...
you loaded a smaller amount of green sample in relation to the red sample.	the peaks for the green sample might appear half as tall as those of the red sample.	To make it easier to view both samples on the same scale, increase the dye scale value of the green sample to make the peaks appear similar.

Changing the Dye Scale of an Electropherogram To change the dye scale of an individual electropherogram:

Step	Action
1	Display the electropherograms with legends.
2	Double-click the dye color indicator next to the sample you want to change. The Choose a Dye Scale dialog box appears. 
3	Enter a new scale in the dialog box and click OK. The dye scale for the individual sample changes, and a vertical line appears beside the dye color indicator to signify that it is modified. Note You can change the dye color in the same way from the Results Control window. The dye is scaled each time you open the Results Display window.

Changing the Dye Scale Preferences

To change the dye scale preferences:

Step	Action
1	<p>Choose Preferences from the Settings menu and choose Results Dye Scales from the submenu.</p> <p>The Preferences dialog box appears with the Results Dye Scales pop-up.</p>  <p>The screenshot shows a window titled "Preferences" with a dropdown menu set to "Results Dye Scales". Inside the window, there are five input fields: "B: 1.0", "Y: 1.0", "R: 1.0", "G: 1.0", and "R: 1.0". Below these fields is a "Reset to Factory Settings" button. At the bottom right of the window are "Cancel" and "OK" buttons.</p>
2	<p>Enter a positive number between 0.1 and 100 for each sample relative to any other sample, and click OK.</p> <p>Note Dye scale values do not automatically revert to default values. Change them back to the defaults before examining results of another run.</p>

Process of Verifying Results

Introduction You can use the electropherogram and tabular displays to verify the results of analysis by checking the GeneScan Analysis Software calculated sizes and peaks.

Note The size calling of the standard and of sample fragments depend on the size calling method you defined in the Analysis Parameters and the accuracy of the defined standard.

Steps to Verifying Size Calculation

To verify size calculations:

Step	Action	See
1	Compare how well multiple size standard electropherograms line up within a Results Display window when aligned by size.	"How to Verify Size Calculations" on page 7-48.
2	View the sizing curve calculated by the GeneScan Analysis Software.	"Size Curve View" on page 3-13.
3	Determine how well the defined size standard matched the size standard run with your sample. Use the Peak Total information in the Sample Info view of the Sample File window.	"Sample Info View" on page 3-8.
4	View the Analysis Log, which provides messages for each analyzed Sample file. If there is a problem or a questionable condition during size calling, a warning message is displayed in the Analysis Log.	"Using the Analysis Log" on page 7-51
5	Use the Raw Data view to display information about the raw data for a sample. Analyzed Sample files contain both raw and analyzed data.	"Raw Data View" on page 3-15.

Step	Action	See
6	<p data-bbox="517 142 864 224">Use EPT Data to troubleshoot problems caused by poor run conditions, such as the:</p> <ul data-bbox="517 240 864 527" style="list-style-type: none"><li data-bbox="517 240 864 300">◆ EP voltage (for ABI PRISM 377).<li data-bbox="517 316 864 341">◆ EP current.<li data-bbox="517 357 864 381">◆ EP power (for the ABI 373).<li data-bbox="517 397 864 457">◆ Laser power (for the ABI PRISM 310).<li data-bbox="517 474 864 527">◆ Gel temperature versus time. <p data-bbox="517 544 864 597">EPT data can be displayed for each Sample file.</p>	"EPT Data View" on page 3-17.

How to Verify Size Calculations

Introduction This section describes the following:

- ◆ Verifying for the GeneScan-350 Standard.
 - ◆ Evaluating for Multiple Internal Size Standards.
-

Verifying for the GeneScan-350 Standard

About the Standard

In the GeneScan-350 standard, the size of the first peak should be approximately 50 bp, the second 75 bp, and so on, assuming that processing started after the 35 bp fragment passed the scan region. If this appears to be correctly measured for your run, measurement of the sample fragments that ran with the standard should also be correct.

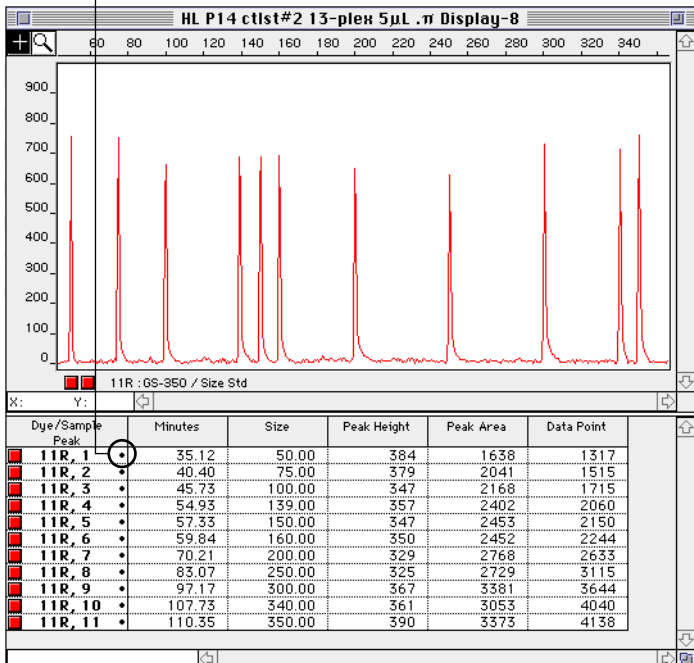
Note For a complete list of fragment sizes, refer to Appendix B, “GeneScan Size Standards.”

Procedure

To verify the size calculation for the GeneScan-350 standard:

Step	Action
1	In the Results Control window, select the Electropherogram and the Tabular Display icons.
2	Click the Clear All button to clear the panels.
3	Click the dye color indicator for the size standard you want to view.
4	Click Display. Note You could also open the Sample File window of the Sample file of interest to verify size calling.

To verify the size calculation for the GeneScan-350 standard: *(continued)*

Step	Action																																																																								
5	<p data-bbox="517 178 1223 235">Click each peak in the electropherogram and check the tabular data to ensure it is the correct size.</p> <p data-bbox="517 251 1169 316">Note The peaks matched to the defined size standard are identified by dots next to the Dye/Sample Peak field.</p> <p data-bbox="517 341 1041 373">Indicates peak matched to defined size standard</p>  <table border="1" data-bbox="517 812 1142 1039"> <thead> <tr> <th>Dye/Sample Peak</th> <th>Minutes</th> <th>Size</th> <th>Peak Height</th> <th>Peak Area</th> <th>Data Point</th> </tr> </thead> <tbody> <tr><td>11R, 1 •</td><td>35.12</td><td>50.00</td><td>384</td><td>1638</td><td>1317</td></tr> <tr><td>11R, 2 •</td><td>40.40</td><td>75.00</td><td>379</td><td>2041</td><td>1515</td></tr> <tr><td>11R, 3 •</td><td>45.73</td><td>100.00</td><td>347</td><td>2168</td><td>1715</td></tr> <tr><td>11R, 4 •</td><td>54.93</td><td>139.00</td><td>357</td><td>2402</td><td>2060</td></tr> <tr><td>11R, 5 •</td><td>57.33</td><td>150.00</td><td>347</td><td>2453</td><td>2150</td></tr> <tr><td>11R, 6 •</td><td>59.84</td><td>160.00</td><td>350</td><td>2452</td><td>2244</td></tr> <tr><td>11R, 7 •</td><td>70.21</td><td>200.00</td><td>329</td><td>2768</td><td>2633</td></tr> <tr><td>11R, 8 •</td><td>85.07</td><td>250.00</td><td>325</td><td>2729</td><td>3115</td></tr> <tr><td>11R, 9 •</td><td>97.17</td><td>300.00</td><td>367</td><td>3381</td><td>3644</td></tr> <tr><td>11R, 10 •</td><td>107.73</td><td>340.00</td><td>361</td><td>3053</td><td>4040</td></tr> <tr><td>11R, 11 •</td><td>110.35</td><td>350.00</td><td>390</td><td>3373</td><td>4138</td></tr> </tbody> </table>	Dye/Sample Peak	Minutes	Size	Peak Height	Peak Area	Data Point	11R, 1 •	35.12	50.00	384	1638	1317	11R, 2 •	40.40	75.00	379	2041	1515	11R, 3 •	45.73	100.00	347	2168	1715	11R, 4 •	54.93	139.00	357	2402	2060	11R, 5 •	57.33	150.00	347	2453	2150	11R, 6 •	59.84	160.00	350	2452	2244	11R, 7 •	70.21	200.00	329	2768	2633	11R, 8 •	85.07	250.00	325	2729	3115	11R, 9 •	97.17	300.00	367	3381	3644	11R, 10 •	107.73	340.00	361	3053	4040	11R, 11 •	110.35	350.00	390	3373	4138
Dye/Sample Peak	Minutes	Size	Peak Height	Peak Area	Data Point																																																																				
11R, 1 •	35.12	50.00	384	1638	1317																																																																				
11R, 2 •	40.40	75.00	379	2041	1515																																																																				
11R, 3 •	45.73	100.00	347	2168	1715																																																																				
11R, 4 •	54.93	139.00	357	2402	2060																																																																				
11R, 5 •	57.33	150.00	347	2453	2150																																																																				
11R, 6 •	59.84	160.00	350	2452	2244																																																																				
11R, 7 •	70.21	200.00	329	2768	2633																																																																				
11R, 8 •	85.07	250.00	325	2729	3115																																																																				
11R, 9 •	97.17	300.00	367	3381	3644																																																																				
11R, 10 •	107.73	340.00	361	3053	4040																																																																				
11R, 11 •	110.35	350.00	390	3373	4138																																																																				
6	<p data-bbox="517 1079 1189 1112">If you find that the sizes were not calculated correctly, you can:</p> <ul style="list-style-type: none"> <li data-bbox="517 1120 893 1153">◆ Redefine the size standard, or <li data-bbox="517 1161 1216 1226">◆ Change the analysis parameters and re-analyze the affected samples (see “Defining Analysis Parameters” on page 5-18). 																																																																								

Evaluating for Multiple Internal Size Standards

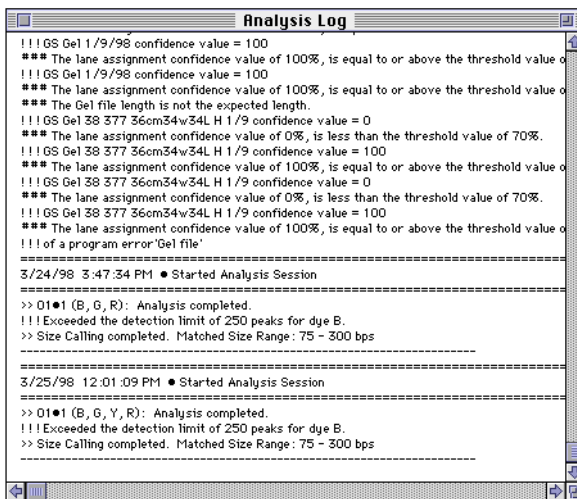
To evaluate size calculations for multiple internal size standards:

Step	Action
1	In the Results Control window, select the Electropherogram button.
2	Click the Clear All button to clear the panels.
3	Click the On radio button to turn on the Quick Tile option.
4	<p>Click the dye color indicators for the size standards you want to view.</p> <p>When the Quick Tile option is on, the GeneScan Analysis Software inserts each in a separate panel.</p> <p>Note Click the header of the appropriate dye/sample column to display all standards in the project that are the same color dye.</p>
5	<p>Click Display.</p> <p>The standards appear in tiled electropherogram displays.</p> <p>For information on setting the tiled electropherogram displays, see “Creating Tiled Electropherogram Displays” on page 7-10.</p>
6	<p>Choose Align by Size (⌘ T) from the View menu if the electropherograms are not already aligned by size.</p> <p>The size standards should line up when aligned by size.</p> <p>Note You can set preferences so that all new displays show data aligned by size.</p> <p>For more information, see “Changing How the Results are Displayed and Printed” on page 7-13.</p>

Using the Analysis Log

What is the Analysis Log The Analysis Log maintains a running record of analysis performed by the GeneScan Analysis Software. If a problem occurs during analysis of a Sample file, the Analysis Log automatically opens and appears in the foreground as an alert.

Displaying the Analysis Log Choose Analysis Log from the Windows menu. The following is an example of the Analysis Log.



What to Evaluate Evaluate the following:

What to evaluate	What not to evaluate
<p>Potential problems that the GeneScan Analysis Software might have had during size calling.</p> <p>It will alert you to any peaks defined in the size standard definition that the software failed to match to the size standard run with your sample.</p>	<p>The GeneScan Analysis Software will not alert you to any consecutive peaks at the end of the definition.</p> <p>This is to avoid logging warnings when your sample was not run long enough to include all the defined size standard peaks.</p> <p>This prevents you from having to create a new standard for shorter runs.</p> <p>The Analysis log will, however, alert you if less than 50% of the defined size standard peaks were not matched, regardless of the peak locations in the definition.</p>

Removing Information from the Analysis Log

To remove information from the Analysis Log:

Step	Action
1	<p>Select the information you want to remove.</p> <p>Note Choose Select All (⌘ A) from the Edit menu to select all the information.</p>
2	<p>Choose Clear from the Edit menu.</p>

Closing the Analysis Log

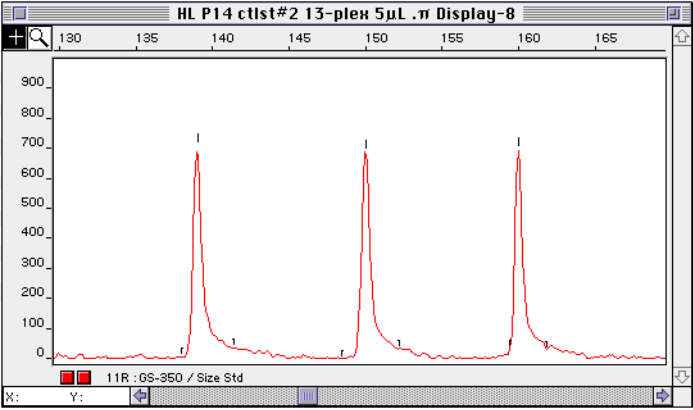
You can either:

- ◆ Click the close box in the upper-left corner.
 - ◆ Choose Close (⌘ W) from the File menu.
-
-

How to Verify Peak Detection

Introduction Use the Show Peak Position command, while the electropherogram and associated tabular data are displayed, to verify results by examining how the GeneScan Analysis Software defined the total area that comprises each peak and the center of the peak.

Verifying Peak Detection To verify peak detection:

Step	Action
1	<p>Choose Show Peak Positions from the View menu.</p> <p>Markers appear that identify the beginning, center, and end of each peak.</p> <p>Note For a better view, choose Zoom In from the View menu, or use the Zoom tool (see “Zooming In and Out” on page 7-31).</p> 
2	<p>Examine the display to ensure that each peak's center, beginning, and end points are correct.</p>
3	<p>Choose Hide Peak Positions from the View menu to suppress the display of the peak markers.</p> <p>Note You can also use the Sample Info view to display information about the peaks detected and matched.</p> <p>For more information, see “Description of Information” on page 3-9.</p>

Saving, Archiving and Copying Files

8

Introduction

In This Section Topics in this chapter include the following:

Topics	See page
Saving GeneScan File	8-2
How to Archive Sample and Gel Files	8-6
Transferring Data to Other Applications	8-7

Saving GeneScan File

Why Save The following table lists why to save projects, Sample files, and gel files. For information on archiving files, see “How to Archive Sample and Gel Files” on page 8-6.

Save	Because	See
GeneScan projects	Protects the links to Sample files and their preferences. Projects contain links to Sample files and preferences regarding display and analysis.	“Saving Projects” on page 8-3.
Sample files	Protects the links to projects and their preferences. Sample files also contain raw data and critical information about the run, settings, and analysis control.	“Saving Sample Files” on page 8-3.
Gel files	Preserves the integrity of the data. Note If you require long-term storage of multiple gel files, saving selected information from the files reduces their size considerably.	“Saving Gel Files” on page 8-4.
Results Displays	Saves the Results Display settings in projects when you have found a display format that suits your needs.	“Saving Results Displays” on page 8-5.

Saving Projects **Note** You do not need to save a Sample file after analysis. The analyzed data is written directly to the Sample file during analysis.

The following table lists the options for saving a project:

If you choose...	Then...						
Save Project (⌘ S)	you can take the following action:						
	<table border="1"><thead><tr><th>If you...</th><th>Then...</th></tr></thead><tbody><tr><td>previously saved the project</td><td>it is automatically saved using the same name.</td></tr><tr><td>had not saved the project</td><td>a dialog box appears. Select a location for the file, enter a name, and click Save.</td></tr></tbody></table>	If you...	Then...	previously saved the project	it is automatically saved using the same name.	had not saved the project	a dialog box appears. Select a location for the file, enter a name, and click Save.
	If you...	Then...					
previously saved the project	it is automatically saved using the same name.						
had not saved the project	a dialog box appears. Select a location for the file, enter a name, and click Save.						
Save Project As	a dialog box appears. Select a location for the file, enter a name, and click Save.						

Saving Sample Choose Save Sample Info (⌘ S) from the File menu.

Files **Note** If you choose Close from the File menu, or click the close box when you have not saved the changes, a dialog box asks if you want to save them.

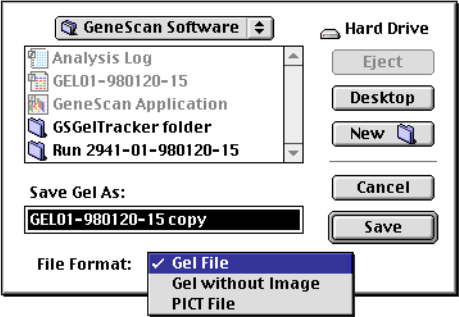
Saving Gel Files **Why Save Only Parts of the Gel File Information**

A gel file can be large, normally 10 MB to 60 MB. You do not need to keep a gel file once the tracking is verified and the corresponding Sample files are created. The following procedure describes how to keep parts of the information, while discarding the rest.

IMPORTANT Do not discard any gel file until you have verified the tracking and taken any required corrective action.

Procedure

To save selected information about a gel file:

Step	Action
1	<p>Select Save As from the File menu.</p> <p>The Save As dialog box appears.</p> 
2	<p>Select a location and type in a name for the file.</p> <p>The default file name is the original file name with “copy” added.</p>

To save selected information about a gel file: *(continued)*

Step	Action								
3	You have the following options from the pop-up menu:								
	<table border="1"><thead><tr><th>File format</th><th>Description</th></tr></thead><tbody><tr><td>Gel file</td><td>Saves the entire gel file into the specified folder. This includes the image, raw data, and other collection information.</td></tr><tr><td>Gel without image</td><td>Saves all of the information in the gel file excluding the image. Storage requirements are about one-third less than the entire gel file. When this gel file is later opened again in the analysis software, the image will be regenerated.</td></tr><tr><td>PICT file</td><td>Creates a PICT file from the gel image. This PICT file can not be opened in the analysis software.</td></tr></tbody></table>	File format	Description	Gel file	Saves the entire gel file into the specified folder. This includes the image, raw data, and other collection information.	Gel without image	Saves all of the information in the gel file excluding the image. Storage requirements are about one-third less than the entire gel file. When this gel file is later opened again in the analysis software, the image will be regenerated.	PICT file	Creates a PICT file from the gel image. This PICT file can not be opened in the analysis software.
	File format	Description							
	Gel file	Saves the entire gel file into the specified folder. This includes the image, raw data, and other collection information.							
Gel without image	Saves all of the information in the gel file excluding the image. Storage requirements are about one-third less than the entire gel file. When this gel file is later opened again in the analysis software, the image will be regenerated.								
PICT file	Creates a PICT file from the gel image. This PICT file can not be opened in the analysis software.								

Saving Results Displays

You can combine electropherograms and tabular data in many ways for display, and the GeneScan Analysis Software allows you to save display combinations and formats for future viewing.

For information on saving a display for future viewing, see “Saving and Renaming the Results Control Format” on page 7-17.

How to Archive Sample and Gel Files

Introduction Since the gel file is so large, you will probably want to back up only specific Sample files. Archive Sample files when you feel confident that the channel selections (tracking) used to generate them was correct.

If you need to save a gel file, you can save only selected information from it, and greatly reduce the file's size (see "Saving Gel Files" on page 8-4).

Archiving Sample and Gel Files

The following table lists how to archive Sample and gel files:

Item	Description	How to archive
Sample files	A sample file is 60 KB to 150 KB in size, depending on the length of the run.	Drag the file icon or the Run folder containing the files to floppy disk icon or to an alternative storage device. A 1.4 MB high-density disks holds about twelve files.
Gel files	Gel files contain the raw data acquired by the Data Collection software and take up a lot of space: 10 MB to 60 MB is typical. After the GeneScan Analysis Software generates Sample files from a correctly-tracked gel file, you do not need to save the gel file.	Make a backup copy of the gel file and remove the file from the hard disk. For best performance, keep as few files on your hard disk as possible. Use magnetic tapes, removable cartridge drives, or optical disks to archive gel files. Gel files are too big for diskettes.

Transferring Data to Other Applications

Other Applications GeneScan files can be read by the following version of the Genotyper® software for further analysis:

Genotyper software version...	Can read...
1.0	Sample files from the GeneScan® Analysis Software
2.0	both Sample files and project files

Both programs can convert GeneScan Analysis Software data into formats used by downstream applications like the GenBase™ software, GenoPedigree™ software, and Linkage, among others.

Cutting and Pasting Tabular Data

To cut and paste tabular data:

Step	Action								
1	Display the tabular data you want to copy.								
2	Select the rows you want to copy by taking the following action:								
	<table border="1"><thead><tr><th>If you want to select...</th><th>Then...</th></tr></thead><tbody><tr><td>all tabular data</td><td>choose Select All (⌘ A) from the File menu.</td></tr><tr><td>several consecutive rows</td><td>shift-click the first and last row in the group you want to select.</td></tr><tr><td>several rows that are not listed next to each other</td><td>⌘-click the rows.</td></tr></tbody></table>	If you want to select...	Then...	all tabular data	choose Select All (⌘ A) from the File menu.	several consecutive rows	shift-click the first and last row in the group you want to select.	several rows that are not listed next to each other	⌘-click the rows.
	If you want to select...	Then...							
	all tabular data	choose Select All (⌘ A) from the File menu.							
several consecutive rows	shift-click the first and last row in the group you want to select.								
several rows that are not listed next to each other	⌘-click the rows.								
3	Choose Copy from the Edit menu.								
4	Open the new application and click where you want to place the information.								
5	Choose Paste from the Edit menu.								

Creating a Text File

To create a text file from tabular data.

Step	Action
1	Display the tabular data.
2	Choose Export Table from the File Menu.
3	Choose a name and file location in the dialog box and click Save.

9

Printing Results

Introduction

In this Chapter Topics in this chapter include the following:

Topics	See page
About Printing	9-2
Automatically Printing Run Results	9-3
Printing the Gel Image	9-4
Printing Selected Sample Files	9-5
Printing Supporting Information	9-6

About Printing

Ways You Can Print

You can print the:

- ◆ Results of your run automatically at the end of the run or interactively, as selected Sample files or display combinations.
 - ◆ Gel file from the Gel File window, and supplementary information from the Sample Info view.
-

If You Get Unexpected Results

You might initially get unexpected results from autoprinting or the Print One command if you switch printers.

The first time you print after changing the printer configuration, do the following:

First	Second
Choose Print Setup from the File menu and click OK.	Choose Print from the File menu and use the standard Print dialog box.

Automatically Printing Run Results

Introduction You can specify that the results are printed in the Data Collection software or in the GeneScan® Analysis Software.

From the Data Collection Software When you choose automatic printing from the Data Collection software, the GeneScan Analysis Software prints a separate page for each designated Sample file, showing electropherograms and tabular data as specified in the Auto-Analysis Defaults (see step 2 on page 4-6).

Choose automatic printing in the Data Collection software as follows:

On this instrument...	Choose...
ABI PRISM® 310	Auto-Print in the Injection list.
ABI PRISM® 377	Auto-Print in the data collection Run window.

For information about setting up your run, refer to your instrument user manual.

From the GeneScan Analysis Software When you analyze a project separately from the Data Collection software, you can print results automatically as the samples are analyzed.

To print the results automatically as the samples are analyzed:

Step	Action
1	In the Analysis Control window, select the samples you wish to analyze. For more information, see “Analyzing a Sample File” on page 3-19.
2	Select the checkbox labeled Print Results in the Analysis Control window.
3	Click the Print Setup button to specify the samples and the format. For more information, see “Specifying the Format for Printed Results” on page 5-9.
4	Click OK. The results are printed after the results are analyzed.

Printing the Gel Image

About Printing the Gel Image

You can print a gel image any time it is displayed.

Note For best results when printing, adjust the gel contrast to increase the intensity of the gel colors. Depending on your printer, the printed image may be less intense than the screen image.

Procedure

To print the gel image, press Command (⌘)-Shift-3 to get a screen capture of everything displayed on the screen.

Printing Selected Sample Files

Introduction You can print selected Sample files using the Results Control window or by choosing the Sample file.

Setting Printing Options To set the printing options, use the Results Display Preferences dialog box (see “Changing How the Results are Displayed and Printed” on page 7-13).

These options affect the electropherogram height and page breaks. Depending on how you set these options, the format that prints may be different from what is on the screen.

For information on printing saved Results Control formats, see “Working with a Previously Saved Displays” on page 7-18.

From the Results Control Window You print selected samples after analysis, regardless of whether you choose automatic printing.

To print results for selected Sample files after analysis:

Step	Action
1	<p>In the Results Control window, select the dye/samples and format you wish to print.</p> <p>Use the same technique as selecting the format and the dye/samples to display the data.</p> <p>For more information, see “Using the Results Control Window” on page 7-7.</p>
2	<p>Take the following action:</p> <ul style="list-style-type: none">◆ Click Print.◆ Choose Print or Print One from the File menu. <p>Note If you choose Print One the Print dialog box does not appear.</p>

From the File Menu

To print a Sample file from the File menu:

Step	Action
1	Choose Open from the File menu. The Open Existing dialog box appears. Note You can also double-click the sample name in the Finder. If the GeneScan Analysis Software is not running, the software starts and opens the Sample file.
2	Click the Sample file icon.
3	In the dialog box that appears, find and select the Sample file that you want to open.
4	Click Open. The Sample File window appears.
5	Select a display mode and choose Print or Print One from the File menu.

Printing Supporting Information

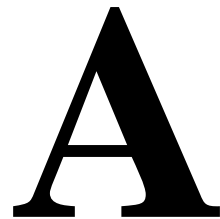
Introduction The following procedure describes how to print supporting information about a Sample file, such as the Sample Info view, or to print the sample sheet.

Note You can use the same procedure for ABI™ 373 and ABI PRISM 377 runs to print the Sample Sheet and an image of the Gel File window.

Procedure To print supporting information:

Step	Action
1	Display the information you want to show.
2	Choose Print (⌘ P) or Print One from the File menu. Note If you choose Print One the Print dialog box does not appear.

Using GeneScan with the ABI 373



Introduction

In This Appendix **Note** This section does not apply if you are processing data collected on ABI™ 373 DNA Sequencer with XL upgrade instruments.

Topics in this appendix include the following:

Topics	See page
Processing Using the ABI 373	A-2
Setting Gel Processing Parameters	A-5
Opening the Collection File	A-7
Completing the Sample Sheet	A-9
Tracking Lanes	A-11
Generating Sample Files	A-12

Processing Using the ABI 373

Introduction This section describes using the ABI 373 data with the GeneScan Analysis Software in the same way as you use ABI PRISM® 377 data. When the GeneScan® 672 Software has been upgraded to work on a Power Macintosh® computer running System 7.5 or higher, this will be true.

**ABI 373 Versus
ABI PRISM 377
Data** The ABI 373 and ABI PRISM 377 run on different computers. If you use the GeneScan® 672 Software to collect data on the ABI 373, the data-collection program runs on a 68K Macintosh® computer. Since the GeneScan Analysis Software 2.0 or later runs on a Power Macintosh computer, transfer the data to a Power Macintosh®-compatible computer before analyzing the data.

**Using the
GeneScan Analysis
Software Before
Version 2.0** Using the GeneScan® 672 Software on data collected using the ABI 373, the software creates a Collection file that contains the raw data. When using previous versions of the GeneScan Analysis Software (before 2.0), you preprocessed this Collection file to create a gel file. Once you were satisfied with the tracking of the gel file, the gel file could be analyzed to create a Results file. The Results file contained all of the analyzed data.

**Using the
GeneScan Analysis
Software Version
2.0 or Later** Using the GeneScan® Analysis Software version 2.0 or later, it was not necessary to preprocess the Collection file. You process the Collection file to create Sample files, then analyze the Sample files. Each Sample file contains information from one lane of the gel and, after analysis, the corresponding analyzed data.

Processing Steps The steps to using the GeneScan Analysis Software with ABI 373 data are as follows:

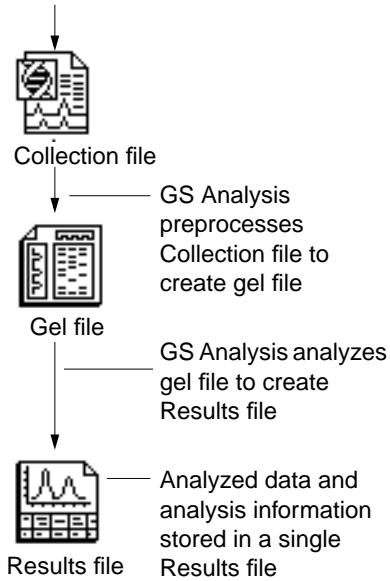
Step	Action	See page
1	Setting Gel Processing Parameters.	A-5
2	Opening the Collection File.	A-7
3	Completing the Sample Sheet.	A-9
4	Tracking lanes.	A-11
5	Generating Sample files.	A-12

Comparison of GeneScan Analysis The following figure is a comparison of using the GeneScan® Analysis Software before version 2.0 and how to use the software with version 2.0 or later. Once you have generated Sample files, use the GeneScan Analysis Software as described in this manual.

Note After analysis, the GeneScan Analysis Software automatically creates a Run folder named “373 SampleFiles <date>”, where <date> is the original date on which the GeneScan® 672 Software file was generated.

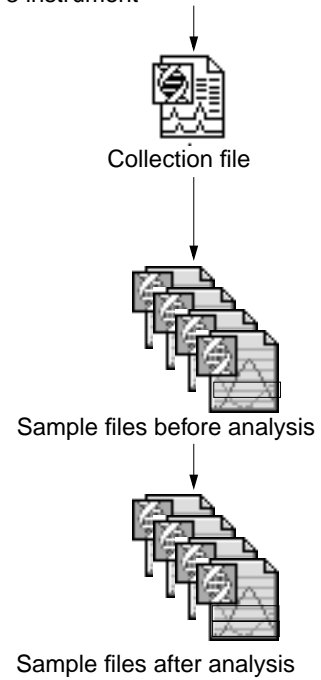
GeneScan Prior to Version 2.0

GeneScan® 672 Software on
ABI 373 instrument



Using GeneScan 2.0 or Later

GeneScan® Analysis Software on
ABI 373 instrument

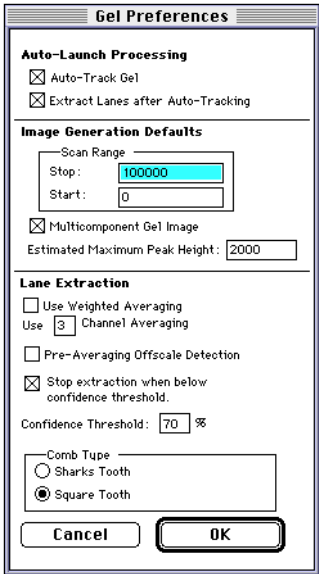


Setting Gel Processing Parameters

Introduction Before you open a Collection file, set the gel processing parameters. The GeneScan® Analysis Software takes the information in the processing parameters into account when it opens the Collection file.

Procedure

To set the gel processing parameters:

Step	Action
1	Double-click the GeneScan Analysis Software icon to start the program.
2	<p>Choose Gel Preferences from the Settings menu.</p> <p>The Gel Preferences dialog box appears.</p> <p>Note The checkboxes under Auto-Launch processing only apply when the Data Collection software automatically starts the GeneScan Analysis Software. The GeneScan® 672 Software does not automatically start GeneScan 2.1.</p> <p>IMPORTANT Before choosing Auto-Track Lane, verify that the comb type is set correctly by choosing Preferences from the Settings menu and Gel Preferences from the submenu.</p> 

To set the gel processing parameters: *(continued)*

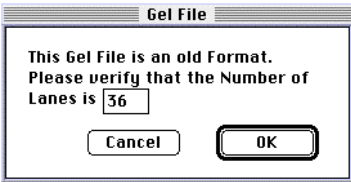
Step	Action								
3	<p>You can take the following action in the Image Generation Defaults section:</p> <table border="1" data-bbox="465 248 1184 708"> <thead> <tr> <th data-bbox="465 248 830 287">To...</th> <th data-bbox="830 248 1184 287">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="465 287 830 435"> start and stop processing data Note All scans outside the specified scan range are ignored. </td> <td data-bbox="830 287 1184 435"> enter the scan numbers in the Scan Range Stop and Start text entry boxes. </td> </tr> <tr> <td data-bbox="465 435 830 505"> compensate for the spectral overlap between dyes </td> <td data-bbox="830 435 1184 505"> select the Multicomponent Gel Image checkbox. </td> </tr> <tr> <td data-bbox="465 505 830 708"> set the upper limit of the scale of the peaks in the Slice view of the Gel window Note This value also affects the brightness of the fragment bands. </td> <td data-bbox="830 505 1184 708"> enter a value in the Estimated Maximum Peak Height text box. </td> </tr> </tbody> </table> <p data-bbox="465 727 1184 784">For more information, see “Image Generation Defaults” on page 2-4.</p> <p data-bbox="465 803 1184 889">Note Setting a range for gel display processing is similar to the same function in the preprocessing parameters of the GeneScan 1.X software.</p>	To...	Then...	start and stop processing data Note All scans outside the specified scan range are ignored.	enter the scan numbers in the Scan Range Stop and Start text entry boxes.	compensate for the spectral overlap between dyes	select the Multicomponent Gel Image checkbox.	set the upper limit of the scale of the peaks in the Slice view of the Gel window Note This value also affects the brightness of the fragment bands.	enter a value in the Estimated Maximum Peak Height text box.
To...	Then...								
start and stop processing data Note All scans outside the specified scan range are ignored.	enter the scan numbers in the Scan Range Stop and Start text entry boxes.								
compensate for the spectral overlap between dyes	select the Multicomponent Gel Image checkbox.								
set the upper limit of the scale of the peaks in the Slice view of the Gel window Note This value also affects the brightness of the fragment bands.	enter a value in the Estimated Maximum Peak Height text box.								
4	<p>You can take the following action in the Lane Extraction section:</p> <table border="1" data-bbox="465 943 1184 1292"> <thead> <tr> <th data-bbox="465 943 830 982">To...</th> <th data-bbox="830 943 1184 982">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="465 982 830 1052"> use a weighted average </td> <td data-bbox="830 982 1184 1052"> select the weighted averaging checkbox. </td> </tr> <tr> <td data-bbox="465 1052 830 1122"> specify the number of channels to be averaged </td> <td data-bbox="830 1052 1184 1122"> enter a value in the Channel averaging text box. </td> </tr> <tr> <td data-bbox="465 1122 830 1292"> stop extraction when the extraction process falls below a set confidence level </td> <td data-bbox="830 1122 1184 1292"> select the Stop extraction when below confidence level checkbox. Enter a value in the Confidence Threshold box. </td> </tr> </tbody> </table> <p data-bbox="465 1312 1184 1341">For more information, see “Lane Extraction” on page 2-7.</p>	To...	Then...	use a weighted average	select the weighted averaging checkbox.	specify the number of channels to be averaged	enter a value in the Channel averaging text box.	stop extraction when the extraction process falls below a set confidence level	select the Stop extraction when below confidence level checkbox. Enter a value in the Confidence Threshold box.
To...	Then...								
use a weighted average	select the weighted averaging checkbox.								
specify the number of channels to be averaged	enter a value in the Channel averaging text box.								
stop extraction when the extraction process falls below a set confidence level	select the Stop extraction when below confidence level checkbox. Enter a value in the Confidence Threshold box.								
5	<p>Click OK.</p>								

Opening the Collection File

Introduction When first opening a Collection file, you usually need to tell the GeneScan Analysis Software how many lanes were collected, and what matrix file should be used to adjust for spectral overlap.

Procedure

Opening the Collection file:

Step	Action
1	Choose Open from the File menu. The dialog box appears.
2	You can either: <ul style="list-style-type: none">◆ Double-click the file name, or◆ Select the desired Collection file and click Open. The gel file verification dialog box appears. 
3	When opening a Collection file, specify to the GeneScan Analysis Software the number of lanes collected by entering the appropriate number in the entry field and clicking OK.
4	Another dialog box appears if the Multicomponent Gel Image checkbox in the Gel Processing Parameters was selected. Identify the matrix file to use for adjusting spectral overlap. For more information, see Chapter 7, "Evaluating Analysis Results." Note If multicomponent was not chosen, but later you decide to regenerate the image with adjustment for spectral overlap, choose Install New Gel Matrix from the Gel menu to attach a Gel matrix (see "Installing New Matrix Information" on page 2-24).

Opening the Collection file: *(continued)*

Step	Action
5	<p data-bbox="471 183 1157 207">Find and select the matrix file you wish to attach and click Open.</p> <p data-bbox="471 232 1177 342">Note This version of the GeneScan Analysis Software can read a matrix generated using GeneScan 1.X. Move the matrix file from the 68K Macintosh® computer to the Power Macintosh® computer you are using for analysis.</p> <p data-bbox="471 367 838 391">The GeneScan Analysis Software:</p> <ul data-bbox="471 407 1089 521" style="list-style-type: none"><li data-bbox="471 407 942 431">◆ Attaches the matrix to the Collection file<li data-bbox="471 448 861 472">◆ Adjusts for spectral overlap, and<li data-bbox="471 488 1089 513">◆ Generates a gel image, which appears on the screen. <p data-bbox="471 537 1049 561">For more information on the gel image, see page 2-27.</p>

Completing the Sample Sheet

Introduction After opening the Collection file, complete the Sample Sheet embedded and track the lanes. This ensures that when the Sample files are generated they contain the correct information.

The Sample Sheet that is part of the Collection file is blank when it is opened. This is because the GeneScan® 672 Software does not automatically embed a copy of the data collection Sample Sheet in the Collection file.

Procedure **IMPORTANT** When tracking and extracting the Collection file, ensure the Sample Sheet associated with the gel file has the checkbox labeled “Used” selected for each sample you want to extract. The GeneScan Analysis Software only extracts the samples with that checkbox selected.

To complete the Sample Sheet:

Step	Action								
1	Open the Sample Sheet by either: <ul style="list-style-type: none"> ◆ Clicking the Sample Sheet button in the upper-left corner of the Gel Window, or by ◆ Choosing Gel Sample Sheet from the Gel menu. 								
2	Enter the following information: <ul style="list-style-type: none"> ◆ Sample file name. ◆ Sample name for each sample. ◆ Sample info. ◆ Comment for each dye/sample. For more information, see “Displaying the Sample Sheet” on page 2-29.								
3	Select checkboxes, as applicable. <table border="1" style="width: 100%; margin-top: 10px;"> <thead> <tr> <th>To...</th> <th>Then select the...</th> </tr> </thead> <tbody> <tr> <td>indicate that the lane contains data</td> <td>Used checkbox.</td> </tr> <tr> <td>automatically analyze Sample files when generating them</td> <td>"A" checkbox.</td> </tr> <tr> <td>automatically print Sample files when generating them</td> <td>"P" checkbox.</td> </tr> </tbody> </table>	To...	Then select the...	indicate that the lane contains data	Used checkbox.	automatically analyze Sample files when generating them	"A" checkbox.	automatically print Sample files when generating them	"P" checkbox.
To...	Then select the...								
indicate that the lane contains data	Used checkbox.								
automatically analyze Sample files when generating them	"A" checkbox.								
automatically print Sample files when generating them	"P" checkbox.								

To complete the Sample Sheet: *(continued)*

Step	Action
4	Identify the dye standard for each sample by clicking in the appropriate dye cells under the standard column. Note Alternately, you can create a stand-alone Sample sheet and attach it to the gel file using the Install New Sample Sheet command from the Gel menu (see "Installing a New Sample Sheet" on page 2-36).
5	Choose Save (⌘-S) or Save As from the File menu. If you click the close box before saving, a dialog box appears with the message whether or not to save the entries.

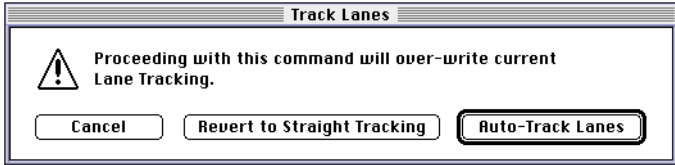
Tracking Lanes

Introduction When opening a Collection file, the GeneScan Analysis Software places an appropriate number of straight tracker lines on the image, based on the number of lanes chosen using the Gel file verification dialog box (see “Opening the Collection File” on page A-7). The software does not automatically track the lanes. You must do so before generating Sample files.

Process The normal process is as follows:

Step	Action
1	Track the lanes.
2	Review and edit the tracking, if necessary.
3	Generate the Sample files. Note However, if you wish, you can automatically extract sample information and create Sample files immediately after tracking.

Procedure To track the lanes and review the tracking before generating Sample files:

Step	Action
1	Choose Track Lanes from the Gel menu. IMPORTANT Before choosing Auto-Track Lane, verify that the comb type is set correctly by choosing Preferences from the Settings menu and Gel Preferences from the submenu. The Track Lanes dialog box appears. 
2	Click Auto-Track Lanes.

Step	Action
3	Review the tracking and edit as necessary. For information on: <ul style="list-style-type: none">◆ Editing tracking, see “Working with Tracker Lines” on page 2-53.◆ Changing the appearance of the window for better visibility, see “How to Adjust the Gel Image” on page 2-18.
4	When you are satisfied with the tracking, choose Save from the File menu.

Generating Sample Files

When to Generate Sample Files

Generate new Sample Files from the lanes if you:

- ◆ Did not automatically generate Sample files after tracking, or you
- ◆ Edited the lane tracking.

For information on generating sample files, see “Extracting Data Without Changing the Current Tracker” on page 2-64.

GeneScan Size Standards

B

Introduction

In This Appendix PE Applied Biosystems provides four different size standards that you can choose from to analyze fragments run on the ABI PRISM® 310, the ABI™ 373, or the ABI PRISM® 377.

See also “About Size Standards” on page 5-29.

Topics in this appendix include the following:

Topics	See page
GeneScan 350 Size Standard	B-2
GeneScan-400HD Size Standard	B-4
GeneScan 500 Size Standard	B-6
GeneScan 1000 Standard	B-8
GeneScan 2500 Standard	B-10

GeneScan 350 Size Standard

What To Use It For GeneScan 350 is a size standard useful for sizing fragments between 35 and 350 base pairs. The native fragments are uniformly spaced to provide accurate size calling.

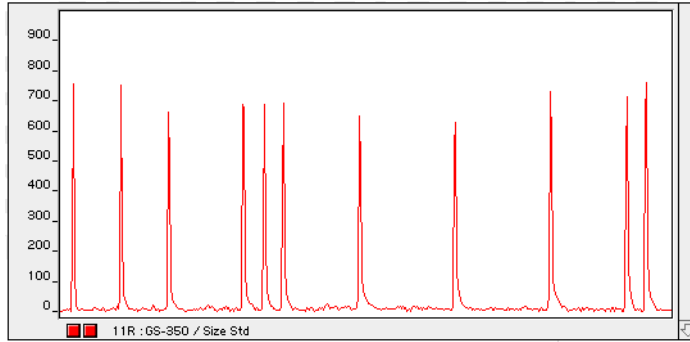
How It Is Prepared GeneScan 350 is prepared by Pst 1 digestion of plasmid DNA, followed by ligation of a TAMRA or ROX-labeled 22-mer oligodeoxynucleotide to the cut ends. A subsequent enzymatic digestion with BstU 1 yields DNA fragments containing a single TAMRA or ROX dye (see “GeneScan 350 Molecular Lengths” below).

GeneScan 350 Molecular Lengths The following table lists the GeneScan 350 Denatured Fragment Molecular Lengths (Nucleotides):

250	139	35
300	150	50
340	160	75
350	200	100

Running Under Denaturing Conditions The following table describes running the GeneScan 350 standard under denaturing conditions:

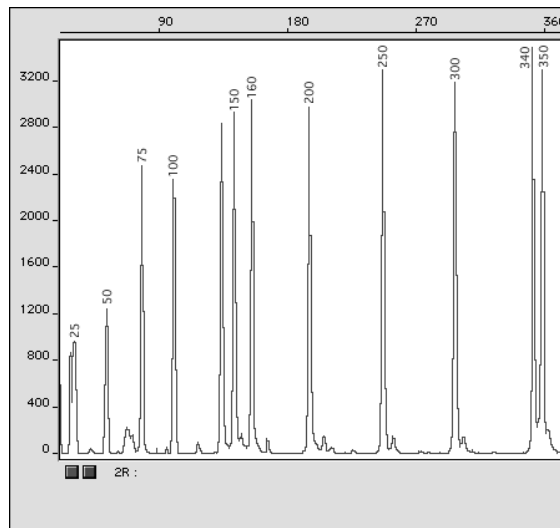
Like	However	Consequently
the GeneScan 2500 and GeneScan 1000 standards, the GeneScan 350 standard is made of double-stranded DNA fragments.	like GeneScan 500, the GeneScan 350 standard has only one labeled strand.	under denaturing conditions, even if the two strands migrate at different rates, only the labeled strand is detected. Because of this, split peaks are avoided that result when two strands move through a denaturing polymer at different rates, see the figure below.



**Double-Stranded
GeneScan 500
Fragments**

The following figure shows the sizes of double-stranded GeneScan 500 fragments. Use these values to size fragments run under native conditions.

IMPORTANT An asterisk (*) for the 250 and 340 base pair peaks denotes peaks resulting from abnormal migration of double strands that did not completely separate under denaturing conditions when analyzed on the ABI PRISM 310. Do not use these peaks to size samples. The peaks show smaller values than the actual size of the fragments. Refer to the *GeneScan Reference Guide, Chemistry Reference for the ABI PRISM 310 Genetic Analyzer* (P/N 4303189 rev A) for further details.



GeneScan-400HD Size Standard

What To Use It For You can use the GeneScan-400HD (High Density) size standard to determine fragment lengths between 50 and 400 base pairs.

Special Uses The high density of marker bands in this standard makes it particularly useful for microsatellite analysis. All fragments have been checked for migration that is true to size under a wide variety of run conditions on all ABI PRISM® instruments. There are no anomalous fragments (*e.g.*, the 250-bp fragment in GeneScan-350 on the ABI PRISM 310 Genetic Analyzer).

Note GeneScan-400HD is the recommended size standard for use with the ABI PRISM Linkage Mapping Set Version 2.

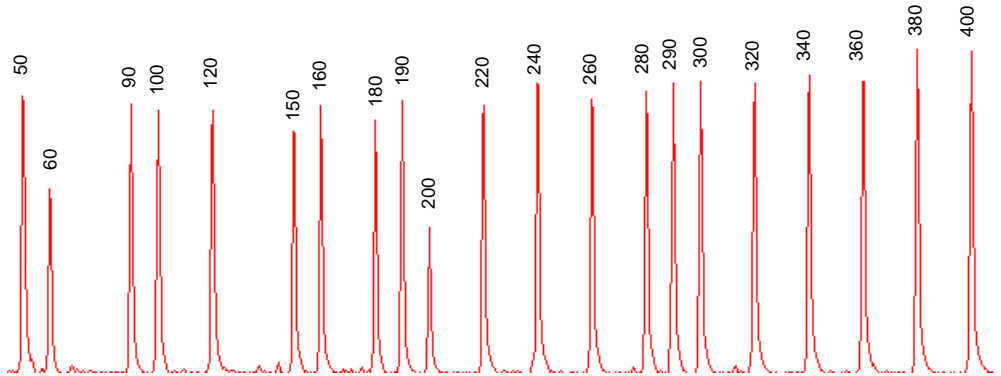
How It Is Prepared All aspects of the preparation of the GeneScan-400HD size standard are proprietary. Each fragment contains a single ROX fluorophore.

Fragment Lengths The following table lists the lengths of the 21 fragments comprising the GeneScan-400HD size standard:

50	160	260	360
60	180	280	380
90	190	290	400
100	200	300	
120	220	320	
150	240	340	

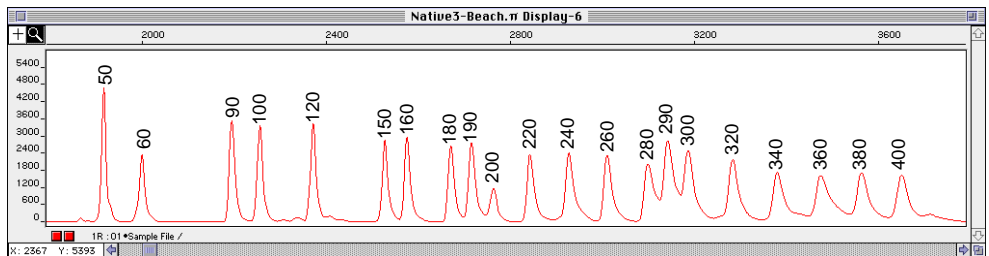
Denaturing Electropherogram

Although the GeneScan-400HD size standard is made of double-stranded DNA fragments, only one of the strands is labeled. Consequently, even if the two strands migrate at different rates under denaturing conditions you will not need to worry about peak splitting. The following figure shows the peak patterns of GeneScan-400HD fragments run under denaturing conditions. Fragments were run using the POP-4 polymer at 60 °C.



Non-denaturing Electropherogram

The following shows the peak patterns of GeneScan-400HD fragments run under non-denaturing conditions. Fragments were run using 3% GeneScan Polymer (GSP) at 30 °C.



GeneScan 500 Size Standard

What To It Use For GeneScan 500 is a size standard useful for sizing fragments between 35 and 500 base pairs. The native fragments are uniformly spaced to provide accurate base calling.

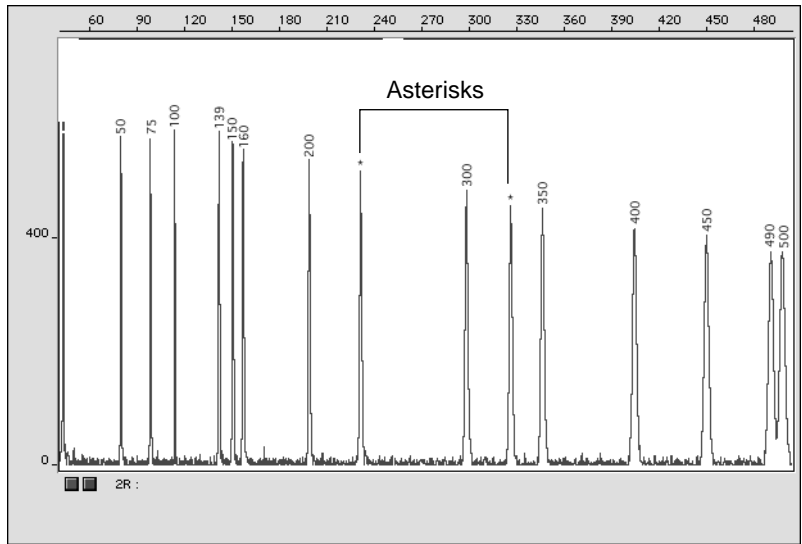
How It Is Prepared GeneScan 500 is prepared by Pst 1 digestion of plasmid DNA, followed by ligation of a TAMRA or ROX-labeled 22-mer oligodeoxynucleotide to the cut ends. A subsequent enzymatic digestion with BstU 1 yields DNA fragments containing a single TAMRA or ROX dye (see “GeneScan 350 Molecular Lengths” below).

GeneScan 500 Molecular Lengths The following table lists the GeneScan 500 Denatured Fragment Molecular Lengths (Nucleotides):

400	250	139	35
450	300	150	50
490	340	160	75
500	350	200	100
400	250	139	35

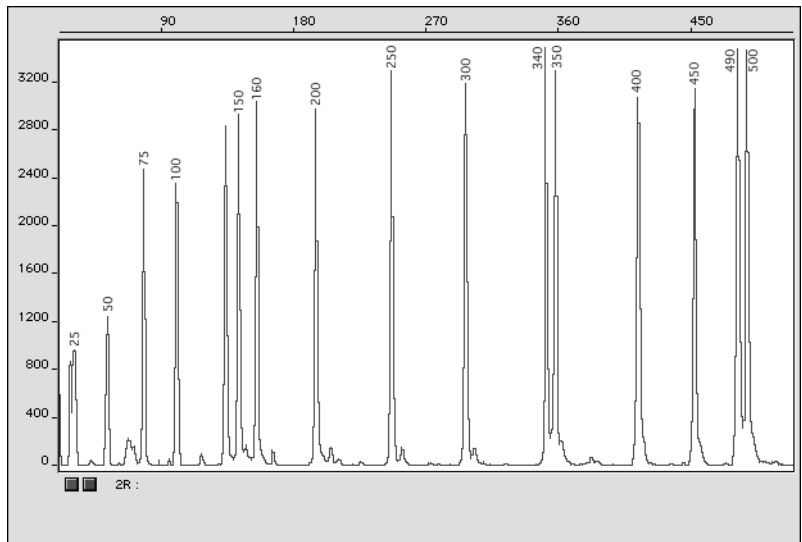
Running Under Denaturing Conditions The following table describes running the GeneScan 500 standard under denaturing conditions:

Like	However	Consequently
the GeneScan 2500 and GeneScan 1000 standards, the GeneScan 500 standard is made of double-stranded DNA fragments.	with GeneScan 500, only one strand of the double-stranded DNA is labeled, whereas the other two standards have labels on both strands.	under denaturing conditions, even if the two strands migrate at different rates, only the one labeled strand is detected. Because of this, split peaks are avoided that result when two strands move through a denaturing polymer at different rates, see the figure below



**Double-Stranded
GeneScan 500
Fragments**

The following figure shows the sizes of double-stranded GeneScan 500 fragments. Use these values to size fragments run under native conditions.



GeneScan 1000 Standard

How It Is Prepared The GeneScan 1000 standard is prepared by digesting pBR322 with the restriction enzyme Alu I, followed by the ligation of a ROX labeled oligonucleotide. This standard has 17 fragments, ranging from 47 to 946 base pairs.

GeneScan 1000 Molecular Lengths Fluorescently labeled native fragments are 18 nucleotides longer than denatured fragments (see the table below). You can use this standard to size fragments in the 100 to 900 base pair range.

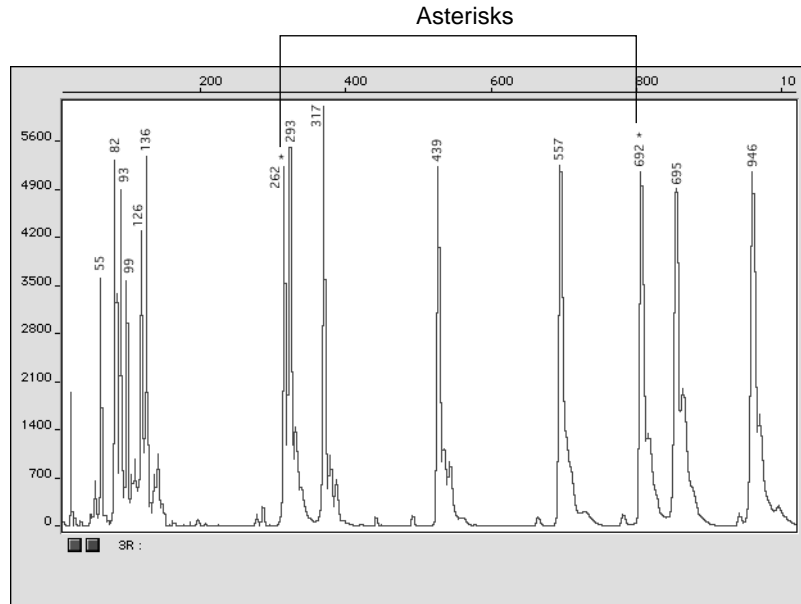
GeneScan 1000 Fragment Molecular Lengths (Base Pairs):

Unlabeled	Native (+36)
928	946
677	695
674	692
539	557
421	439
299	317
275	293
244	262
118	136
108	126
81	99
75	93
56	85
64	82
37	55
33	51
29	47

Running Under Native Conditions

The following diagram shows the peak pattern of fragments run under native conditions on an ABI PRISM 310 using 2.5% GeneScan Polymer Solution in a 30 cm Ld capillary.

IMPORTANT An asterisk (*) for the 262 and 692 base pair peaks denotes peaks resulting from abnormal migration of double strands that did not completely separate under denaturing conditions when analyzed on the ABI PRISM 310. Do not use these peaks to size samples. The peaks show smaller values than the actual size of the fragments. Refer to the *GeneScan Reference Guide, Chemistry Reference for the ABI PRISM 310 Genetic Analyzer* (P/N 4303189 rev A) for further details.



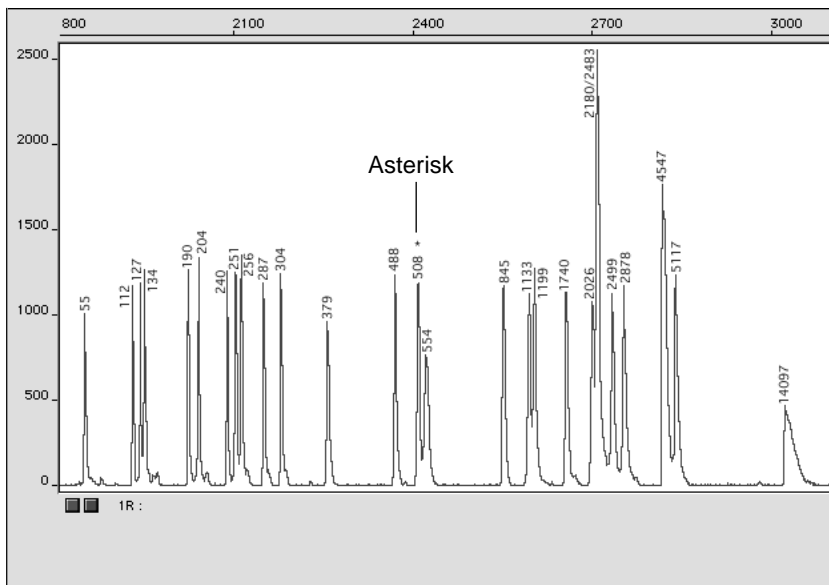
GeneScan 2500 Standard

How It Is Prepared The GeneScan 2500 standard is made from lambda phage DNA restriction digested with Pst I, followed by ligation of either a TAMRA or ROX oligonucleotide. It has 28 fragments, ranging from 55 to 14,097 base pairs (bp).

What To Use It For You can use the GeneScan 2500 standard for native applications to size fragments in the 100 to 5,000 base pair range.

Running Under Native Conditions The following figure shows the peak pattern of fragments run under native conditions on the ABI PRISM 310 using 2.5% GeneScan Polymer Solution in a 30 cm Ld capillary.

IMPORTANT An asterisk (*) for the 508 base pair peaks denotes peaks resulting from abnormal migration of double strands that did not completely separate under denaturing conditions when analyzed on the ABI PRISM 310. Do not use these peaks to size samples. The peaks show smaller values than the actual size of the fragments. Refer to the *GeneScan Reference Guide, Chemistry Reference for the ABI PRISM 310 Genetic Analyzer* (P/N 4303189 rev A) for further details.



**GeneScan 2500
Molecular Lengths**

Fluorescently labeled native fragments are 18 nucleotides longer than denatured fragments (see table below).

GeneScan 2500 Molecular Lengths (bp)

Denatured	Native (+18)
14079	14097
5099	5117
4771	4789
4529	4547
2860	2878
2481	2499
2465	2483
2162	2180
2008	2026
1722	1740
1181	1199
1115	1133
827	845
536	554
490	508
470	488
361	379
286	304
269	287
238	256
233	251
222	240
186	204
172	190
116	134
109	127
94	112
37	55

Size Calling Methods

C

Introduction

In This Appendix Topics in this appendix includes the following:

Topic	See
Least Square Method	C-2
Cubic Spline Interpolation Method	C-4
Local Southern Method	C-5
Global Southern Method	C-7

Least Square Method

About This Method Both Least Squares methods (2nd Order and 3rd Order) use regression analysis to build a best-fit size calling curve. This curve compensates for any fragments that may run anomalously. As a result, this method normally results in the least amount of error for all the fragments, including the size standards and the samples.

Depending on whether you choose the 2nd or 3rd Order Least Squares Method in the Analysis Parameters dialog box, the resulting size curve is either a quadratic or a cubic function. The software uses the known standard fragments and the associated scan number positions to produce a sizing curve based on Multiple Linear Regression.

Advantages In the figures Figure C-1 and Figure C-2, you can see that in nearly all instances the mobility of an individual DNA fragment is coincident with the best curve fit of the entire data set. Stated differently, the mobility of most DNA fragments is strictly length-dependent. This method automatically compensates for fragments that run anomalously.

GeneScan® Analysis Software calculates a best-fit least squares curve for all samples, regardless of the size calling method you choose. The curve is black in the Standard Sizing Curve window.

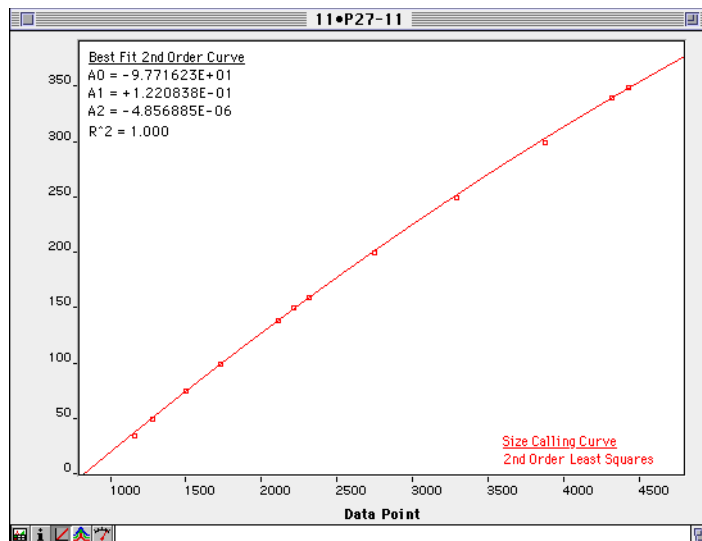


Figure C-1 2nd Order Least Squares size calling curve

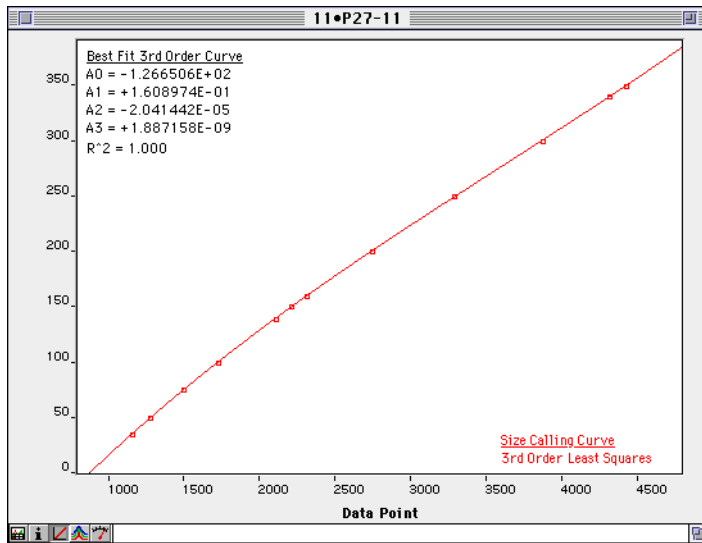
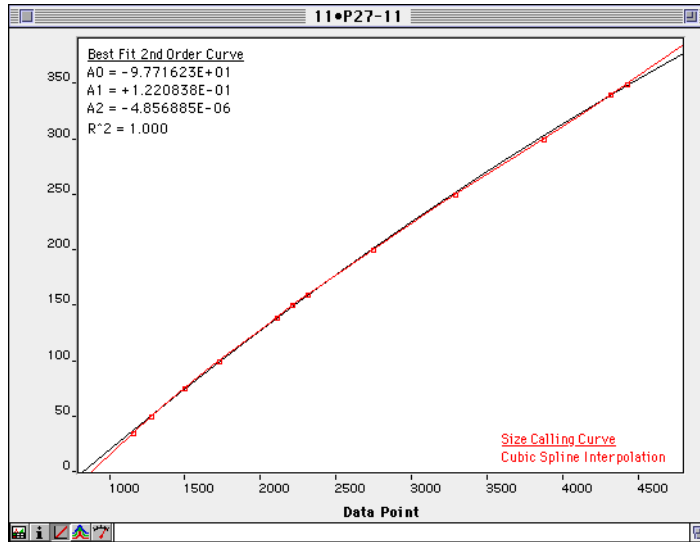


Figure C-2 3rd Order Least Squares size calling curve

Cubic Spline Interpolation Method

About This Method By definition, the Cubic Spline method forces the sizing curve through all the known points of the selected GeneScan size standard. Although this produces exact results for the values of the standards themselves, it does not compensate for standard fragments that may run anomalously.



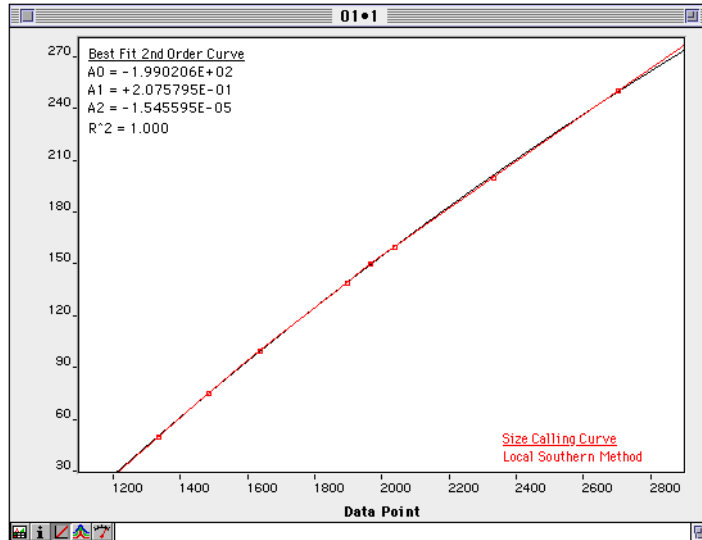
Possible Local Sizing Inaccuracy Mobility of any DNA fragment can be affected by its sequence, and by secondary and tertiary structure formation. If any internal size standard fragment has anomalous mobility, the Cubic Spline method may exhibit local sizing inaccuracy.

For example: Assume that a standard fragment is close in molecular length to an unknown sample fragment. Assume further that the standard fragment runs anomalously. The Cubic Spline method assigns the official value to this standard fragment, even though it may be slightly incorrect. The size of the unknown fragment is then likely to be calculated incorrectly as well.

Note This method does not determine the amount of sizing accuracy error.

Local Southern Method

About This Method The Local Southern method determines the sizes of fragments by using the reciprocal relationship between fragment length and mobility, as described by E. M. Southern (1979).



The Equation The following table describes how the equation works:

Equation	Description
$L = [c/(m-m_0)] + L_0$	Attempts to describe the reciprocal relationship between the mobility, m , and the length, L_0 , of the standard fragments.

How This Method Works

This method, which is similar to the Cubic Spline method, uses the four fragments closest in size to the unknown fragment to determine a best fit line value. Using this method, only the region of the size ladder near the fragment of unknown length is analyzed.

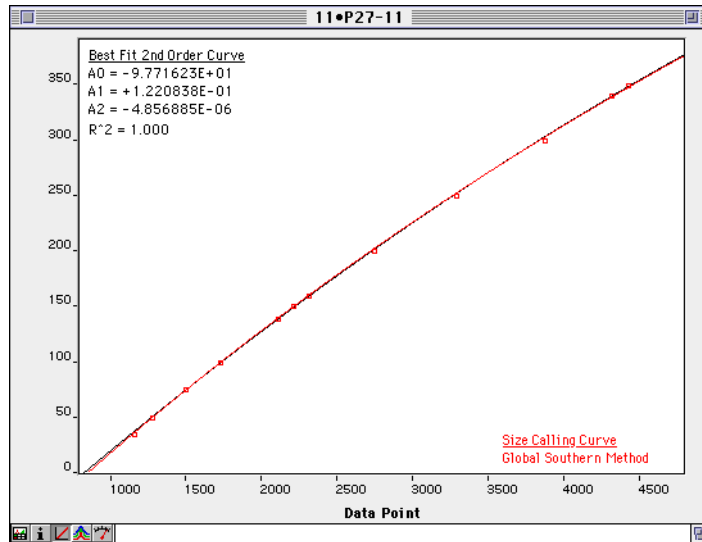
Note Size estimates may be off if any of the standard fragments run anomalously.

The following table lists how the Local Southern method works:

Step	Action
1	The fitting constants of the curve are calculated for each group of three neighboring points on the standard. A separate curve is created for each set of three points.
2	A curve is then created by using three standard points (two points below and one point above the fragment) and a fragment size is determined.
3	Another curve is created by looking at an additional set of three points (one point below and two points above the fragment) and another value is assigned.
4	The two size values are averaged to determine the unknown fragment length.

Global Southern Method

About This Method This method is similar to the Least Squares method in that it compensates for standard fragments that may run anomalously. The method creates a best-fit line through all the available points, and then uses values found on that line to calculate the fragment values.



The Equations The following table describes the equations work:

Equation	Description
$L = [c/(m-m_0)] + L_0$	Attempts to describe the reciprocal relationship between the mobility, m , and the length, L_0 , of the standard fragments.
$\sum_i (L_i - (c/(m_i - m_0) + L_0))^2$	The fitting constants L_0 , m_0 and c are calculated by a least squares fit to minimize the following quantity.

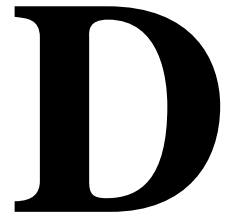
How This Method Works

All points in the standard are weighted equally and the curve is not constrained to go through any specific point. The software can analyze a large range of fragment sizes with this method.

DNA fragments that are...	Are sized using...
not bracketed within the size standard curve	a second order least squares curve extrapolation.
bracketed within the size standard curve	the method that was chosen.

For best results, use a standard that brackets all the fragments of interest.

Troubleshooting the GeneScan Software



Introduction

In This Appendix The tables in this section present information about problems you might experience with your GeneScan® Analysis Software runs, and suggest possible causes and corrections.

For information on troubleshooting the BioLIMS™ database, see Appendix F, “Troubleshooting the BioLIMS Database.”

Topics in this appendix include the following:

Topics	See page
Troubleshooting Projects and Results	D-2
Troubleshooting Gel Data	D-5
Troubleshooting Genotyping Results	D-8
GeneScan Error Messages	D-9

Troubleshooting Projects and Results

Table Description The following table lists the problem, probable cause, and correction for troubleshooting projects and results.

Troubleshooting projects and results:

Problem	Probable Cause	Correction
File name is dimmed in Size Standard or Parameters column in Analysis Control window.	The file has been moved from the folder in which it was located when it was first selected.	<ol style="list-style-type: none"> Move the file back to its original location. Reset preferences to specify the new folder location. Create or select a new file.
Peaks appear on display but the GeneScan Analysis Software does not detect them (cannot select them in electropherogram display).	<ul style="list-style-type: none"> ◆ Peak Amplitude Threshold set too high. ◆ Minimum Peak Half Width set too high. ◆ Electrophoresis run too quickly resulting in poor resolution. 	<ol style="list-style-type: none"> Adjust minimum peak height to include smallest peaks desired and re-analyze. Reduce minimum peak half width setting and re-analyze. Repeat electrophoresis at reduced power. <p>For more information, see “Peak Detection Parameter Options” on page 5-20.</p>
At the position of one strong peak additional colors appear underneath the peak.	<ul style="list-style-type: none"> ◆ Off-scale data not multicomponented correctly. ◆ Poor / incorrect matrix. 	<ol style="list-style-type: none"> Repeat electrophoresis; load less sample. Attach a new gel matrix and regenerate Sample files, or assign a new matrix to the Sample file and re-analyze.

Troubleshooting projects and results: *(continued)*

Problem	Probable Cause	Correction
Peaks appearing in a dye color that should not be present.	Bleed-through from other colors because of off-scale data.	Repeat electrophoresis; load less sample.
A sample run on two different gels does not give the same molecular weights.	<ul style="list-style-type: none"> ◆ Size standard peaks called incorrectly by software. ◆ Different sizing method used for each gel. ◆ Inaccuracies associated with off-scale data do not allow for identification of top of true peak. 	<ul style="list-style-type: none"> a. Check size standard and sizing curve; Re-analyze the Sample file with a different size standard or create a new one. b. Make sure same sizing method has been used for both gels. c. Repeat electrophoresis; load less sample.
Using GS-350 standard, size peaks >350 bp cannot be sized.	<p>Sizing requires at least one standard fragment larger than peak to be sized.</p> <p>For more information, see Appendix B, "GeneScan Size Standards."</p>	Repeat electrophoresis using GS-500.
Peak centers seem to be incorrect in electropherogram.	<ul style="list-style-type: none"> ◆ Resolution of the gel might be inadequate (ABI™ 373 or ABI PRISM® 377). ◆ Signal-to-noise ratio might be too low. 	Repeat electrophoresis at lower power or with longer gel.

Troubleshooting projects and results: *(continued)*

Problem	Probable Cause	Correction
Software cannot display the sizing curve for a sample.	<ul style="list-style-type: none"> ◆ Sample's in-lane size standard does not match defined size standard. ◆ Sample file was not size called. 	<ul style="list-style-type: none"> ◆ Re-analyze the Sample file with a different size standard, or create a new one. <p>In the Analysis Control window:</p> <ul style="list-style-type: none"> d. ⌘-click dye/sample that represents Size Standard for the Sample file. e. Choose a size standard definition file and re-analyze.
Peaks disappear in the electropherogram.	Included the primer peak in the analysis.	

Troubleshooting Gel Data

Table Description The following table lists the problem, probable cause, and correction for troubleshooting gel data.

Troubleshooting gel data:

Problem	Probable Cause	Correction
At the position of one strong peak additional colors appear underneath the peak.	<ul style="list-style-type: none"> ◆ Off-scale data not multicomponented correctly. ◆ Poor / incorrect matrix. ◆ Gel Image not multicomponented. 	<ol style="list-style-type: none"> a. Repeat electrophoresis; load less sample. b. Attach a new gel matrix, regenerate the gel image. c. Regenerate the gel image with multicomponenting selected.
Peaks appearing in a dye color that should not be present.	Bleed-through from other colors because of off-scale data.	Repeat electrophoresis; load less sample.
TAMRA-labeled size standard appears yellow on the gel display.	Collected using: <ul style="list-style-type: none"> ◆ Filter set A (ABI 373), or ◆ Virtual filter A (ABI PRISM 377). 	Repeat electrophoresis with: <ul style="list-style-type: none"> ◆ Filter set B (ABI 373), or ◆ Virtual filter C (ABI PRISM 377).
TET-labeled products not seen on gel display.	Collected using: <ul style="list-style-type: none"> ◆ Filter set A (ABI 373), or ◆ Virtual filter A (ABI PRISM 377). 	Repeat electrophoresis with: <ul style="list-style-type: none"> ◆ Filter set B (ABI 373), or ◆ Virtual filter C (ABI PRISM 377).

Troubleshooting gel data: (continued)

Problem	Probable Cause	Correction
Signal showing up in neighboring lanes.	Leaking wells of gel.	<ul style="list-style-type: none"> ◆ Consider using a square tooth comb instead of a shark tooth comb. ◆ If using 96-lanes, then rerun gel using protocol in the <i>ABI PRISM 377 DNA Sequencer 96-Lane Upgrade User's Manual</i> (P/N 4305423).
	Signal intensity very high and signal is being detected in neighboring lanes due to closeness of spacing.	Move tracker lane position from center of band to edge of the band away from strong signal and extract as usual.
		Use 1 or 2 lane averaging to extract lanes.
HEX-labeled products appear green on gel display.	Collected using: <ul style="list-style-type: none"> ◆ Filter set A (ABI 373), or ◆ Virtual filter A (ABI PRISM 377). 	Repeat electrophoresis with: <ul style="list-style-type: none"> ◆ Filter set B (ABI 373), or ◆ Virtual filter C (ABI PRISM 377).
Collection time was sufficient, but only a small portion of gel displayed.	<ul style="list-style-type: none"> ◆ Gel Image Processing preferences did not include enough scans to display entire gel. ◆ Electrophoresis power too low. 	<ol style="list-style-type: none"> a. Regenerate gel image with new scan range. b. Adjust to correct settings; repeat electrophoresis.

Troubleshooting gel data: *(continued)*

Problem	Probable Cause	Correction
Improper tracking results.	Bad matrix.	Attach new matrix.
	Sample Sheet not filled out properly.	Fill out Sample Sheet properly.
	Comb types set improperly.	a. Fix and type in gel preferences. b. Retrack gel.
	Peak height or red signal too low.	Rerun gel with more size standard.

Troubleshooting Genotyping Results

Table Description The following table lists the problem, probable cause, and correction for troubleshooting the Genotyper® software results.

Problem	Probable Cause	Correction
Allele peaks seen in correct molecular weight range, with additional peaks seen outside this range.	<ul style="list-style-type: none"> ◆ Bleed-through from other colors because of off-scale data. ◆ Primers not fully optimized. 	<ul style="list-style-type: none"> a. Repeat electrophoresis; load/inject less sample. b. Check optimization.
With allele peaks of high intensity, the GeneScan Analysis Software calls many small peaks.	<ul style="list-style-type: none"> ◆ Background above minimum peak height. ◆ Too much PCR product loaded. 	<ul style="list-style-type: none"> a. Adjust minimum peak height; re-analyze. b. Repeat electrophoresis; load/inject less sample. <p>For more information, see "Peak Detection Parameter Options" on page 5-20.</p>
A homozygous individual shows a dip at the top of an allele peak which may be called as two separate peaks.	Truncated single peak because of off-scale data can appear as two peaks.	Repeat electrophoresis; load/inject less sample.
Warning message: "Could not complete 'Run Macro' command because the labeled peak could not be found."	The first allele peak for one or more loci in the allelic ladder is lower than the preset minimum peak height specification in the categories list.	<ul style="list-style-type: none"> a. Adjust minimum peak height; re-analyze. b. Repeat electrophoresis; load/inject less sample. <p>For more information, see "Peak Detection Parameter Options" on page 5-20.</p>

GeneScan Error Messages

Introduction This section includes three tables:

- ◆ GeneScan Analysis Software Crashes with BioLIMS.
- ◆ Analysis Log Error Messages.
- ◆ Error Messages When Defining Size Standards.

GeneScan Analysis Software Crashes with BioLIMS

The following table lists the error message you might encounter if the application crashes with BioLIMS installed:

Observation	Possible Cause	Recommended Action
GeneScan Analysis Software (with BioLIMS) crashes on launch or gives an error message, 'Could not open " " '.	Insufficient memory to load the Oracle® or Sybase® libraries.	Ensure that there is at least 2 MB of free memory in addition to the preferred memory requirements set on the GeneScan Analysis Software Info box. To access the Info box, select the GeneScan application and choose Get Info from the File menu.
GeneScan Analysis Software (with BioLIMS) crashes when attempting to connect to a Sybase database.	SybaseConfig control panel or libtcp extension file is missing or disabled.	Use the Extension Manager to check that these files are present and turned on. If either file is missing, then reinstall the BioLIMS Client software.

Analysis Log Error Messages

The following table lists the error messages you might encounter in the GeneScan Analysis Software Log.

Analysis Log error messages:

Analysis Log Error Message	Comment/Correction	Refer To
The Analysis Range does not include enough data points. Check your Analysis Parameters.	Make sure the Analysis Range in your Analysis Parameters contains at least 250 data points.	“Defining Analysis Parameters” on page 5-18.
The Range of Data Points to analyze is too large. Check your Analysis Parameters.	Specify a smaller range in the Analysis Parameters.	“Defining Analysis Parameters” on page 5-18.
The Analysis Parameters could not be accessed. Check your Analysis Parameters Setting.	Make sure the Analysis Parameters file specified in the Analysis Control window is valid and accessible.	“Defining Analysis Parameters” on page 5-18.
The Sample File does not contain a valid matrix. Assign a new matrix to the Sample File and try again.	Assign a new matrix to the Sample file and re-analyze.	“Installing a New Matrix File” on page 3-20.

Analysis Log error messages: *(continued)*

Analysis Log Error Message	Comment/Correction	Refer To
<p>The Sample File's Matrix was not found.</p> <p>Assign a new matrix to the Sample File and try again.</p>	<p>Assign a new matrix to the Sample file and re-analyze.</p> <p>Choose Preferences from the Settings menu and Folder Locations from the submenu.</p> <p>Check the location of the matrix folder.</p> <p>For more information, see "Specifying File Locations" on page 5-42.</p>	<p>"Installing a New Matrix File" on page 3-20.</p>

Error Messages When Defining Size Standards

The following table list the error messages you might encounter while defining size standards.

Error messages when defining size standards:

Error Message	Comment/Correction	Refer To
<p>The affected Sample File is not available.</p> <p>Locate the Sample File and try again.</p>	<p>If the Sample file name is dim in the Analysis Control window, the GeneScan Analysis Software has not located the sample file.</p> <p>You can instruct the program to search for the sample file.</p>	<p>"Finding Missing Sample Files" on page 4-14.</p>
<p>A Dye Standard is not selected for the affected Sample File.</p> <p>Select a Dye Standard and try again.</p>	<p>Select the dye/sample that represents the standard by ⌘-clicking the appropriate dye/sample field.</p>	<p>Callout 3 on page 5-3.</p>

Error messages when defining size standards: *(continued)*

Error Message	Comment/Correction	Refer To
<p>The affected Sample File does not have a valid Analysis Parameters Selection.</p> <p>Select new Analysis Parameters and try again.</p>	<p>Select either:</p> <ul style="list-style-type: none"> ◆ The default program parameters (<Analysis Parameter >), or ◆ A valid analysis parameters file in the Analysis Control window. 	<p>Callout 5 on page 5-3.</p>
<p>No peaks were found within the Analysis Range.</p> <p>Check your Analysis Parameters.</p>	<ul style="list-style-type: none"> ◆ Make sure the Peak Amplitude Threshold setting allows for detection of the peaks in your sample. ◆ If peaks in your data are narrow, make sure the Minimum Peak Half Width is a small number. 	<p>“Defining Analysis Parameters” on page 5-18.</p>

Using the BioLIMS Database



Introduction

In This Appendix This appendix describes how to access the BioLIMS™ database, how to set the preferences, and how to open or process a fragment that is located in the BioLIMS database.

For information on troubleshooting the BioLIMS database, see Appendix F, “Troubleshooting the BioLIMS Database.”

Topics in this appendix include the following:

Topic	See page
About the GeneScan Analysis Software and the BioLIMS Database	E-2
Configuring the BioLIMS Database Server	E-4
Switching Between Sample File and BioLIMS Mode	E-10
How to Access the BioLIMS Database	E-12
About Server Names	E-17
Using the Collection Browser Window	E-20

About the GeneScan Analysis Software and the BioLIMS Database

What is the BioLIMS Database? The GeneScan® Analysis Software 3.1 introduced support for selecting, viewing, editing, and processing fragments from the BioLIMS database.

The BioLIMS database provides a relational database for fragments created by ABI PRISM® GeneScan Analysis Software. This database accommodates multiple users and versions while preserving the original data.

Accessing the BioLIMS Database The GeneScan Analysis Software user interface has changed to accommodate accessing the BioLIMS database.

Once you have accessed the BioLIMS database, you can display the Collection Browser window by choosing the Open command from the File menu.

Before Using the BioLIMS Database Before you can connect to the BioLIMS database you must have installed the BioLIMS 2.0 software. It is recommended that you install the BioLIMS software first, then install the GeneScan Analysis Software into the BioLIMS 2.0 software folder.

Note This organization allows you to install the Sequencing Analysis software at a later time.

For more information, see “Installing the Software” on page 1-12.

Modes in GeneScan Analysis Software If you are using the GeneScan Analysis Software with the BioLIMS database, you can use the software in either of two modes:

Mode	Descriptions
Sample File mode	In Sample File mode, fragment data extracted from gel files is written out to individual sample files.
BioLIMS mode	In BioLIMS mode, fragment data extracted from gel files is written to a BioLIMS database that resides on a Oracle® Server database or Sybase SQL Server™. Sample files that came from the ABI PRISM 310 Genetic Analyzer can be uploaded to the BioLIMS database using the Sample 2DB Software.

Comparing Modes The following table is a comparison of BioLIMS and Sample File modes.

Feature	BioLIMS mode	Sample File mode
data extracted from a gel file is written to	the BioLIMS database	sample files
fragment data is opened for viewing and for analysis from	the BioLIMS database	sample files

Configuring the BioLIMS Database Server

Sybase or Oracle? This section gives instructions on how to configure the client computer (that runs the GeneScan Analysis Software) for database access.

The BioLIMS database resides on either a Sybase SQL Server or an Oracle Server.

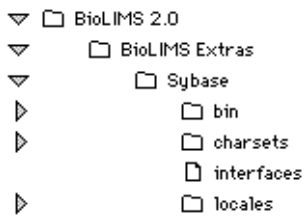
To configure for the...	See...
Sybase SQL Server	"Configuring for Sybase SQL Server Connection" on page E-4.
Oracle Server	"Configuring for Oracle Server Connection" on page E-7.

Configuring for Sybase SQL Server Connection

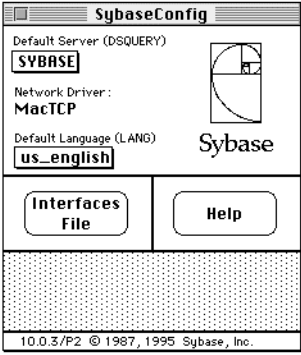
Follow the steps below to configure a Macintosh® computer for connection to the Sybase SQL Server.

IMPORTANT Any time the name, port number, IP address, or host and domain name of the BioLIMS database server is changed, you need to repeat this procedure.

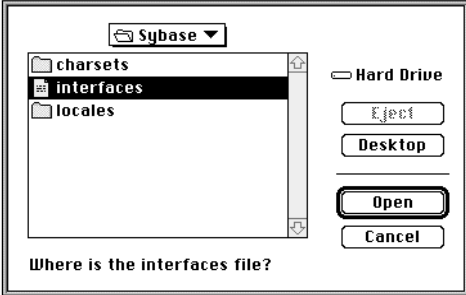
To configure for Sybase SQL Server connection:

Step	Action
1	Find the interfaces file in the Sybase® folder in the BioLIMS Extras folder.  <pre>▼ [] BioLIMS 2.0 ▼ [] BioLIMS Extras ▼ [] Sybase ▶ [] bin ▶ [] charsets ▶ [] interfaces ▶ [] locales</pre>
2	Open the file with SimpleText or a similar text editing application.

To configure for Sybase SQL Server connection: *(continued)*

Step	Action
3	<p>Find the lines:</p> <pre>SYBASE query MacTCP mac_ether neuron.apldbio.com 2500</pre> <p>and edit them:</p> <ul style="list-style-type: none"> ◆ Replace SYBASE by an alias name for the database server (see “About Server Names” on page E-17). ◆ Replace neuron.apldbio.com with the IP address or host and domain name of the server machine. ◆ 2500 is the default port number for the Sybase database. If necessary, replace 2500 with the port number recommended by your BioLIMS database administrator. <p>You can find this information in the interfaces file on the Sybase Server, or your BioLIMS database administrator can provide the information.</p>
4	<p>If you have access to more than one server, duplicate the two lines and edit them for the other server(s).</p> <p>For example, for two servers, one called SYBASE and one called SERVER2, the interfaces file might look like this:</p> <pre>SYBASE query MacTCP mac_ether neuron.apldbio.com 2500 SERVER2 query MacTCP mac_ether 192.135.191.128 2025</pre>
5	<p>Save and close the interfaces file.</p>
6	<p>Open the SybaseConfig control panel. This control panel is found in the Control Panels folder in the System folder.</p> 

To configure for Sybase SQL Server connection: *(continued)*

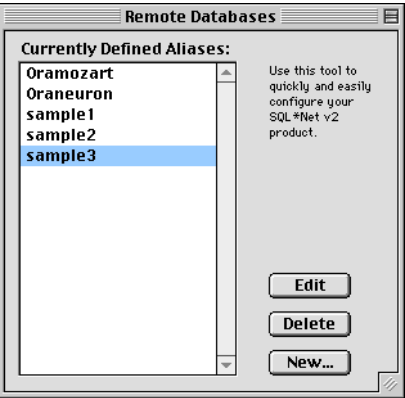
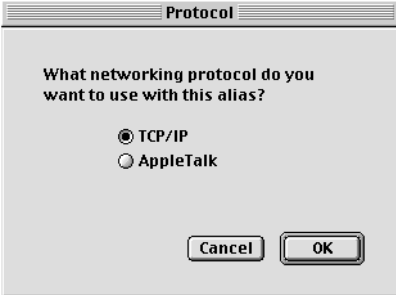
Step	Action
7	<p>The first time the SybaseConfig control panel is opened, a file browser opens automatically.</p> <p>If a file browser does not open immediately, click the Interfaces Files button to open one.</p> <div data-bbox="475 347 939 639" style="border: 1px solid black; padding: 5px;"><p>The screenshot shows a file browser window with a title bar that says 'Sybase'. Inside the window, there is a list of folders: 'Charsets', 'interfaces', and 'locales'. The 'interfaces' folder is selected and highlighted. To the right of the list, there are several buttons: 'Hard Drive', 'Eject', 'Desktop', 'Open', and 'Cancel'. The 'Open' button is highlighted with a thick border. Below the list, there is a text prompt: 'Where is the interfaces file?'.</p></div>
8	Use the file browser to locate and open the interfaces file edited in the steps above.
9	Set the Default Language pop-up menu to us_english.
10	Close the SybaseConfig control panel.

Configuring for Oracle Server Connection

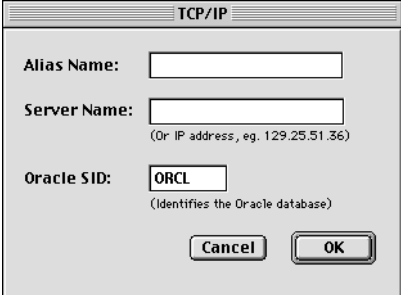
Follow the steps below to configure your Macintosh computer for connection to the Oracle Server.

IMPORTANT Any time you change the BioLIMS database server name or its IP address or host (and domain name), you need to repeat this procedure.


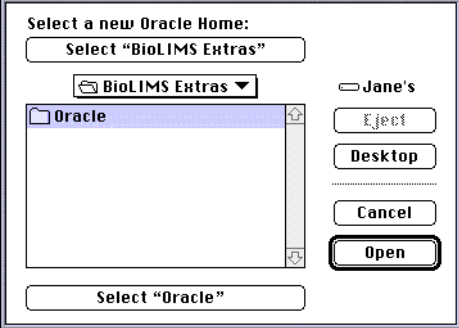
To configure for Oracle Server connection:

Step	Action
1	<p>Use the program Easy Config to configure for the Oracle Server connection.</p> <p>At installation, Easy Config is placed into the BioLIMS 2.0:BioLIMS Extras:Oracle:Applications:Networking folder.</p> <p>Open the Easy Config program.</p> <p>The Remote Databases window appears.</p> 
2	<p>Click New.</p> <p>The Protocol dialog box appears.</p> 

To configure for Oracle Server connection: *(continued)*

Step	Action
3	<p>Select TCP/IP and click OK.</p> <p>The TCP/IP dialog box appears.</p> 
4	<p>Enter text in the fields as follows:</p> <ul style="list-style-type: none"> ◆ Alias Name: Enter an alias name for the database server. (See “About Server Names” on page E-17.) ◆ Server Name: Enter the server name. This may be an IP address or host (and domain name) of the server machine. <p>Note This field does not scroll horizontally for display even though it accepts characters typed past the end of the field. If the server name is longer than 20 characters, you may want to enter the end characters first and go back or just use the IP address.</p> <ul style="list-style-type: none"> ◆ Oracle SID: Enter the value of the ORACLE_SID environment variable. <p>You can find this information in the tnsnames.ora file on the Oracle Server, or your BioLIMS database administrator can provide you with the information.</p>
5	Click OK to close the TCP/IP dialog box.
6	From the File menu, choose Save Configuration.
7	From the File menu, choose Quit to exit the Easy Config program.

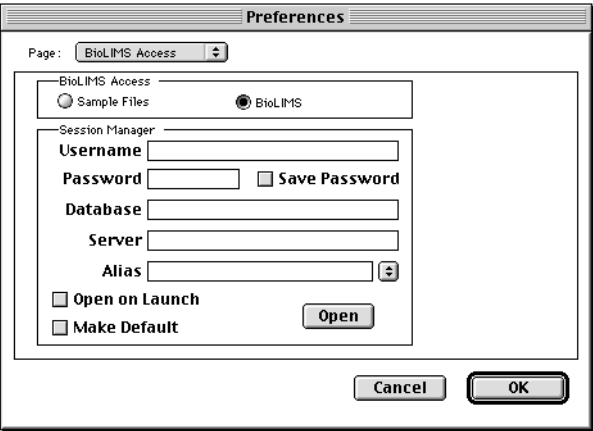
To configure for Oracle Server connection: *(continued)*

Step	Action
8	<p>Find the application Set Oracle Home.</p>  <p>Set Oracle Home</p> <p>The application is contained in your BioLIMS 2.0:BioLIMS Extras:Oracle:Applications folder.</p> <ul style="list-style-type: none"> ▼ <input type="checkbox"/> BioLIMS 2.0 ▼ <input type="checkbox"/> BioLIMS Extras ▼ <input type="checkbox"/> Oracle ▼ <input type="checkbox"/> Applications <ul style="list-style-type: none"> ◆ Set Oracle Home ◆ SQL *Plus ▶ <input type="checkbox"/> Libraries ▶ <input type="checkbox"/> Network
9	<p>Open the application Set Oracle Home.</p> 
10	<p>Use the file browser to find the Oracle folder in the BioLIMS Extras folder.</p> <p>Press Select Oracle button.</p>

Switching Between Sample File and BioLIMS Mode

Introduction If the BioLIMS database is installed on your system, then the BioLIMS Access command is available from the Settings menu when you start the GeneScan Analysis Software.

Switching Modes To switch modes in the GeneScan Analysis Software:

Step	Action									
1	<p>Choose Preferences from the Settings menu and BioLIMS Access from the submenu.</p> <p>The Preferences dialog box appears.</p> 									
2	<p>To switch modes, take the following action:</p> <table border="1" data-bbox="465 1076 1188 1346"> <thead> <tr> <th data-bbox="465 1076 709 1117">To change to...</th> <th data-bbox="709 1076 951 1117">Then...</th> <th data-bbox="951 1076 1188 1117">Result</th> </tr> </thead> <tbody> <tr> <td data-bbox="465 1117 709 1214">Sample File mode</td> <td data-bbox="709 1117 951 1214">click the Sample Files radio button.</td> <td data-bbox="951 1117 1188 1214">The Session Manager text is greyed out.</td> </tr> <tr> <td data-bbox="465 1214 709 1346">BioLIMS mode</td> <td data-bbox="709 1214 951 1346">click the BioLIMS radio button.</td> <td data-bbox="951 1214 1188 1346">The fields in the Session Manager subpanel are activated.</td> </tr> </tbody> </table>	To change to...	Then...	Result	Sample File mode	click the Sample Files radio button.	The Session Manager text is greyed out.	BioLIMS mode	click the BioLIMS radio button.	The fields in the Session Manager subpanel are activated.
To change to...	Then...	Result								
Sample File mode	click the Sample Files radio button.	The Session Manager text is greyed out.								
BioLIMS mode	click the BioLIMS radio button.	The fields in the Session Manager subpanel are activated.								

Step	Action						
3	<p data-bbox="521 139 874 167">You can take the following action:</p> <table border="1" data-bbox="517 207 1239 386"><thead><tr><th data-bbox="517 207 659 245">Click...</th><th data-bbox="659 207 1239 245">To...</th></tr></thead><tbody><tr><td data-bbox="517 245 659 318">OK</td><td data-bbox="659 245 1239 318">affect the changes to the Preferences dialog box and to close it.</td></tr><tr><td data-bbox="517 318 659 386">Cancel</td><td data-bbox="659 318 1239 386">close the Preferences dialog box without changing the mode.</td></tr></tbody></table>	Click...	To...	OK	affect the changes to the Preferences dialog box and to close it.	Cancel	close the Preferences dialog box without changing the mode.
Click...	To...						
OK	affect the changes to the Preferences dialog box and to close it.						
Cancel	close the Preferences dialog box without changing the mode.						

How to Access the BioLIMS Database

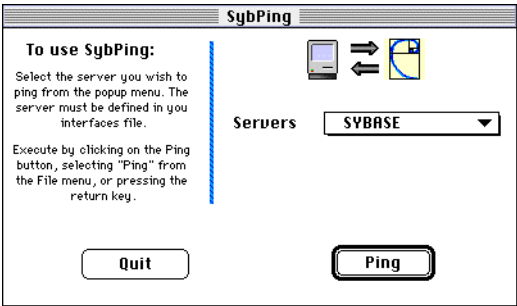
Introduction The following procedure describes how to access the BioLIMS database by completing the Preferences dialog box.

Before you can work with the GeneScan Analysis Software software, you must establish a connection to the BioLIMS database. This connection is made through the BioLIMS Access page of the Preferences dialog box.

Before Accessing a Sybase Database

IMPORTANT If you try to connect to a Sybase database through the GeneScan Analysis Software when the network is down, the computer may hang. If this happens, you have to restart the Macintosh computer. Force the computer to restart by pressing the key combination, Command–Control–Power key.

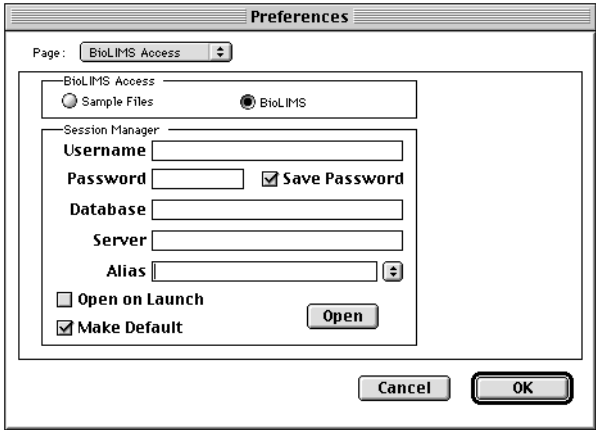
Before accessing a Sybase database, you need to ping the database using SybPing to ensure that the database connection is working.

Step	Action
1	<p>Open the SybPing program.</p> <p>At installation this program is placed in the BioLIMS 2.0:BioLIMS Extras:Sybase:bin folder.</p>  <p>The screenshot shows a Macintosh-style window titled 'SybPing'. On the left, there is text: 'To use SybPing: Select the server you wish to ping from the popup menu. The server must be defined in your interfaces file. Execute by clicking on the Ping button, selecting "Ping" from the File menu, or pressing the return key.' Below this text are two buttons: 'Quit' and 'Ping'. On the right side of the window, there is a 'Servers' label and a dropdown menu currently showing 'SYBASE'. Above the dropdown menu are two icons: a computer icon and a globe icon, with arrows pointing between them.</p>
2	<p>Select the name of the database Server that you want to use.</p> <p>These Server names correspond to those listed in the interfaces file.</p>
3	<p>Click Ping.</p>


Step	Action						
4	<table border="1"> <thead> <tr> <th>If the connection is...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>working</td> <td>the Ping window reports that the "Ping was successful."</td> </tr> <tr> <td>not working</td> <td> <ul style="list-style-type: none"> ◆ The Ping window reports that the "Ping Failed" or, ◆ The Ping window gives no response. <p>Refer to Appendix F, "Troubleshooting the BioLIMS Database."</p> </td> </tr> </tbody> </table>	If the connection is...	Then...	working	the Ping window reports that the "Ping was successful."	not working	<ul style="list-style-type: none"> ◆ The Ping window reports that the "Ping Failed" or, ◆ The Ping window gives no response. <p>Refer to Appendix F, "Troubleshooting the BioLIMS Database."</p>
	If the connection is...	Then...					
	working	the Ping window reports that the "Ping was successful."					
not working	<ul style="list-style-type: none"> ◆ The Ping window reports that the "Ping Failed" or, ◆ The Ping window gives no response. <p>Refer to Appendix F, "Troubleshooting the BioLIMS Database."</p>						
5	Quit from the SybPing application.						

Accessing the BioLIMS Database

To access the BioLIMS database:

Step	Action
1	<p>Choose Preferences from the Edit menu and BioLIMS Access from the submenu.</p> <p>The Preferences dialog box appears.</p> 

To access the BioLIMS database: *(continued)*

Step	Action
2	<p>In the BioLIMS Access section, click the BioLIMS button. The fields in the Session Manager subpanel are activated.</p> <p>Note The following alert appears if there are unsaved sample files and fragments open when you switch to BioLIMS mode. When you click OK, the alert is dismissed and the Preferences dialog box appears with the Sample Files button selected.</p> <div data-bbox="475 404 944 574" style="border: 1px solid black; padding: 5px;"><p>The Sample Data Access mode cannot be changed while editing Sample/Project Files. Please make sure all open Sample/Project Files have been saved.</p><p style="text-align: right;"><input type="button" value="OK"/></p></div>
3	<p>In the Session Manager section in the Preferences dialog box, enter:</p> <ul style="list-style-type: none">◆ Your user name on the server.◆ Your password for your server account.◆ The name of the database on the server (You may have access to more than one database on the server.)◆ The server name. The server name is contained in the interfaces file (Sybase) or in the tnsnames.ora file (Oracle). <p>IMPORTANT All these text boxes are case sensitive.</p>
4	<p>Click the check box labeled Save Password if you want to:</p> <ul style="list-style-type: none">◆ Save your password so that you do not have to enter it every time you open the connection.◆ Run AppleScripts that don't contain password information.◆ Automatically analyze the data after the Data Collection software has collected the data. <p><input checked="" type="checkbox"/> Save Password</p>

To access the BioLIMS database: *(continued)*

Step	Action						
5	<p>If you want the database to open automatically when you launch the GeneScan Analysis Software, click the check box labeled Open on Launch.</p> <p><input checked="" type="checkbox"/> Open on Launch</p> <p>Note You must also click the check box labeled Save Password if you want the database to open automatically in order to automatically analyze data.</p>						
6	<p>If you intend to use more than one database or user account, enter an alias for this BioLIMS session information. You can use an alias to connect to the database if no database connection is open.</p> <p>Once you enter an alias, you can take the following action:</p> <table border="1" data-bbox="517 641 1239 792"> <thead> <tr> <th data-bbox="517 641 705 680">If you click...</th> <th data-bbox="705 641 1239 680">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 680 705 719">Open</td> <td data-bbox="705 680 1239 719">you connect to the database.</td> </tr> <tr> <td data-bbox="517 719 705 792">OK</td> <td data-bbox="705 719 1239 792">the changes are saved and you are connected to the database if a channel is open.</td> </tr> </tbody> </table>	If you click...	Then...	Open	you connect to the database.	OK	the changes are saved and you are connected to the database if a channel is open.
If you click...	Then...						
Open	you connect to the database.						
OK	the changes are saved and you are connected to the database if a channel is open.						
7	Use the pop-up menu to add, change, or remove aliases.						
8	<p>If you have more than one alias, click the Make Default check box to choose which one appears when you first open the Edit Session dialog box.</p> <p><input checked="" type="checkbox"/> Make Default</p> <p>Note The default alias is the database that opens if you choose to automatically analyze data.</p> <p>Note If both the Make Default and the Save Password boxes are checked, no dialog box will appear when a connection to the server is requested. Since all the information required of the user has been saved, the software will connect to the database automatically.</p>						

To access the BioLIMS database: *(continued)*

Step	Action						
9	Take one of the following actions:						
	<table border="1"><thead><tr><th data-bbox="467 253 704 289">If the login was...</th><th data-bbox="704 253 1170 289">Then...</th></tr></thead><tbody><tr><td data-bbox="467 289 704 565">successful</td><td data-bbox="704 289 1170 565">a. Choose Open from the File menu. The Open Existing dialog box appears. b. Click the Sample icon and the Collection Browser appears. For more information, see "Using the Collection Browser Window" on page E-20.</td></tr><tr><td data-bbox="467 565 704 889">unsuccessful</td><td data-bbox="704 565 1170 889">an alert dialog box appears. Check that:<ul style="list-style-type: none">◆ All the login information was entered correctly and in the correct case.◆ Your interfaces file is correctly configured (page E-4).◆ If the connection is still not open, consult Appendix F, "Troubleshooting the BioLIMS Database."</td></tr></tbody></table>	If the login was...	Then...	successful	a. Choose Open from the File menu. The Open Existing dialog box appears. b. Click the Sample icon and the Collection Browser appears. For more information, see "Using the Collection Browser Window" on page E-20.	unsuccessful	an alert dialog box appears. Check that: <ul style="list-style-type: none">◆ All the login information was entered correctly and in the correct case.◆ Your interfaces file is correctly configured (page E-4).◆ If the connection is still not open, consult Appendix F, "Troubleshooting the BioLIMS Database."
	If the login was...	Then...					
successful	a. Choose Open from the File menu. The Open Existing dialog box appears. b. Click the Sample icon and the Collection Browser appears. For more information, see "Using the Collection Browser Window" on page E-20.						
unsuccessful	an alert dialog box appears. Check that: <ul style="list-style-type: none">◆ All the login information was entered correctly and in the correct case.◆ Your interfaces file is correctly configured (page E-4).◆ If the connection is still not open, consult Appendix F, "Troubleshooting the BioLIMS Database."						

About Server Names

Sybase or Oracle? The BioLIMS Session Manager decides whether you are connected to Sybase SQL Server or to an Oracle Server database by looking at the name in the Server field in the Session Manager dialog box.

How Names are Recognized The table below summarizes how names are recognized.

If the Session Manager sees a Server name...	It assumes a...	Example
All in uppercase letters	Sybase SQL Server database connection.	MOZART
Suffixed by ":s" or ":S"	Sybase SQL Server database connection.	Offenbach:S
Containing any lowercase letters	Oracle Server database.	Oramozart
Suffixed by ":o" or ":O"	Oracle Server database.	SIBELIUS:O

Sybase SQL Server Examples

Example 1

If the interfaces file contains this:

```
MOZART
query MacTCP mac_ether mozart.apldbio.com 2500
```

MOZART is recognized as a Sybase SQL Server because the server name is in all uppercase letters.

The Session Manager would look like this:

Username

Password **Save Password**

Database

Server

Example 2

If the interfaces file contains this:

```
Offenbach
query MacTCP mac_ether mozart.apldbio.com 2500
```

The Session Manager would look like this:

Username	<input type="text" value="jane"/>
Password	<input type="password" value="••••••"/> <input type="checkbox"/> Save Password
Database	<input type="text" value="biolims2"/>
Server	<input type="text" value="Offenbach:S"/>

In order for Offenbach to be recognized as a Sybase SQL Server, the name in the Server field is suffixed with "S".

Oracle Server Examples

Example 1

If the tnsnames.ora file contains this:

```
Oramozart=(DESCRIPTION=
            (ADDRESS=
              (PROTOCOL=TCP)(host=mozart)(port=1521))
            (CONNECT_DATA=(SID=WG733))
          )
```

Oramozart is recognized as an Oracle Server because the server name begins with "O".

The Session Manager would look like this:

Username	<input type="text" value="jane"/>
Password	<input type="password" value="••••••"/> <input type="checkbox"/> Save Password
Database	<input type="text" value="biolims2"/>
Server	<input type="text" value="Oramozart"/>

Example 2

If the tnsnames.ora file contains this:

```
SIBELIUS = (DESCRIPTION=
            (ADDRESS=
              (PROTOCOL=TCP) (host=SIBELIUS) (port=1521))
            (CONNECT_DATA=(SID=WG733))
          )
)
```

The Session Manager would look like this:

Username	<input type="text" value="jane"/>
Password	<input type="password" value="••••••"/> <input type="checkbox"/> Save Password
Database	<input type="text" value="biolims2"/>
Server	<input type="text" value="SIBELIUS:O"/>

In order for SIBELIUS to be recognized as a Oracle Server, the name in the Server field is suffixed with “:O”.

Using the Collection Browser Window

Applications that Use the Collection Browser Window

Collection Browser window is common to the following BioLIMS-aware applications.

- ◆ AutoAssembler™ DNA Sequence Assembly Software.
- ◆ GeneScan® Analysis Software.
- ◆ Factura™ Feature Identification Software.
- ◆ Sample 2DB Software.
- ◆ Sequencing Analysis software.

Ways to Search the Database

Using the Collection Browser window from within GeneScan Analysis Software, you can search the BioLIMS database for specific collections and fragments.

The following table lists ways you can search:

Search by...	See page
up to 5 collection-specific criteria	E-24
up to 14 fragment-specific criteria	E-25

In This Section

This section includes the following topics.

For this topic	See page
Displaying the Collection Browser Window	E-21
Collection Browser Window Example	E-22
Parts of the Collection Browser Window	E-22
Collection Search Criteria	E-24
Fragment Search Criteria	E-25
Searching the BioLIMS Database	E-28

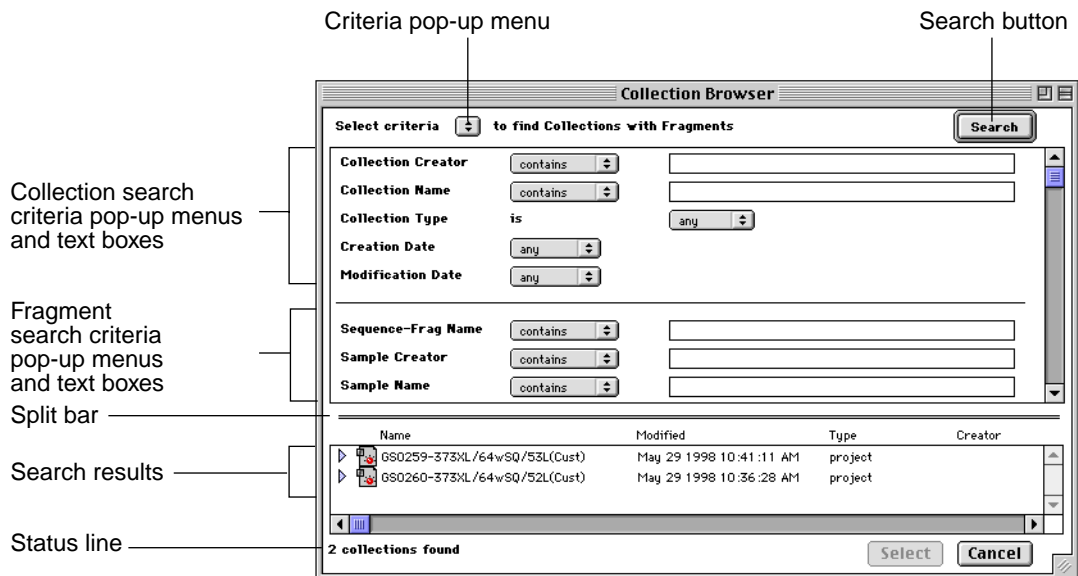
Displaying the Collection Browser Window

There are two ways to display the Collection Browser window:

If you want to...	Then...	Result
open a fragment to view or edit	<ol style="list-style-type: none">Choose Open from the File menu. The Open Existing dialog box opens.Click the Sample icon.	The Collection Browser window appears. For more information, see "Collection Browser Window Example" on page E-22.
create a new project by adding sample files	<ol style="list-style-type: none">Choose New from File menu. The Create New dialog box appears.Click the Project icon. An untitled Analysis Control window appears.Choose Add Samples Files from the Project menu.	

Collection Browser Window Example

The following is an example of the Collection Browser window.



Parts of the Collection Browser Window

The table below describes the parts of the Collection Browser window that were labeled in the figure above.

Item	Description
Criteria pop-up menu	Use this pop-up menu to specify the search criteria visible on the Collection Browser window. Note If you only intend to use a subset of criteria, setting only that subset visible helps to reduce clutter in the window. The search results are the same whether a criterion is invisible or blank and visible.
Search button	Click this button to query the BioLIMS database. Note You can also press the Return key to begin a search.

Item	Description
Collection search criteria pop-up menus and text boxes	<p>Use these pop-up menus and text boxes to define the collection criteria of the search.</p> <p>IMPORTANT Only those fragments that match each and every criterion you specify are returned. That is, search criteria are combined using the logical AND operation.</p> <p>For more information, see “Collection Search Criteria” on page E-24.</p>
Fragment search criteria pop-up menu and text boxes	<p>Use these pop-up menus and text boxes to define the fragment criteria of the search.</p> <p>IMPORTANT A collection is returned if one or more of the fragments contained in it fulfill all of the specified fragment criteria.</p> <p>Note Only fragments meeting search criteria will be displayed in the Collection Browser window.</p> <p>For more information, see “Fragment Search Criteria” on page E-25.</p>
Split bar	<p>Drag this bar to alter the relative amount of space allocated to the top and bottom portions of the Collection Browser window.</p>
Search results	<p>After a successful query, found sample files are listed in this area as Name, Modification date, type, and Creator.</p>
Status line	<p>Search results, error messages, and other important information is reported here.</p> <p>For example, the Status Line lists how many collections were returned in a search.</p>

Collection Search Criteria

The table below shows the collection search criteria. The collections returned by the Collection Browser window must match all of the collection criteria and contain at least one fragment that matches all of the fragment criteria.

Allowed Collection Search Criteria:

Criterion	Pop-up Menu Choices	Allowed Text	Description
Collection Creator	◆ is ◆ starts with ◆ ends with ◆ contains	up to 255 characters	Name of the creator/owner of the collection.
Collection Name	◆ is ◆ starts with ◆ ends with ◆ contains	up to 31 characters	Name of the collection.
Collection Type	◆ any ◆ run ◆ project ◆ other	NA	Collection type. Default is any menu item.
Creation Date	◆ any ◆ is ◆ before ◆ after ◆ between	date — set with arrow buttons using the format mm/dd/yy	Date the collection was created.
Modification Date	◆ any ◆ is ◆ before ◆ after ◆ between	date — set with arrow buttons using the format mm/dd/yy	Date the collection was last modified.

Fragment Search Criteria

The table below shows the fragment search criteria. The collections returned by the Collection Browser window must contain at least one fragment that matches all of the specified fragment criteria.

Fragment Search Criteria

Criterion	Pop-up Menu Choices	Allowed Text	Description
Sequence-Frag Name	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 31 characters including letters, numbers, and punctuation Cannot use colons (:).	Name of the fragment. This is the file name entered in the Sample Sheet.
Sample Creator	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 255 characters including letters, numbers, and punctuation	Name of the person responsible for the run.
Sample Name	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 255 characters including letters, numbers, and punctuation	Sample name from the Sample Sheet.
Instrument Name	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 255 characters including letters, numbers, and punctuation	Set in the General Settings Preferences of the Data Collection software.
Instrumentation	<ul style="list-style-type: none"> ◆ any ◆ gel ◆ capillary 	NA	Whether the sample was run on a gel or capillary instrument.
Start Collect Date	<ul style="list-style-type: none"> ◆ any ◆ is ◆ before ◆ after ◆ between 	date— set with arrow buttons using format mm/dd/yy	Date data collection began.

Fragment Search Criteria *(continued)*

Criterion	Pop-up Menu Choices	Allowed Text	Description
End Collect Date	<ul style="list-style-type: none"> ◆ any ◆ is ◆ before ◆ after ◆ between 	date — set with arrow buttons using format mm/dd/yy	Date data collection ended.
Gel Path	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 255 characters including letters, numbers, and punctuation	The full path name to the original gel file, e.g., Hard Disk:Data: GelRuns:L28t.
Sample Info	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 255 characters including letters, numbers, and punctuation	Sample information from the Sample Sheet.
Sample Comment	<ul style="list-style-type: none"> ◆ is ◆ starts with ◆ ends with ◆ contains 	up to 255 characters including letters, numbers, and punctuation	Comment from the Sample Sheet.
Size Data	<ul style="list-style-type: none"> ◆ is present ◆ is not present ◆ does not apply 	NA	is present means that one or more dyes contain sizing information.
			is not present means none of the dye sample contain sizing information.

Fragment Search Criteria *(continued)*

Criterion	Pop-up Menu Choices	Allowed Text	Description
Size Calling	<ul style="list-style-type: none"> ◆ done ◆ not done ◆ does not apply 	NA	done means sample file has completed size calling indicated by a size curve.
			not done indicated by a missing size curve.
% Matched Peaks	<ul style="list-style-type: none"> ◆ any ◆ equal to ◆ less than ◆ greater than ◆ between 	0—100	Percentage based on size standard matched peaks divided by size standard defined peaks.
Offscale Data	<ul style="list-style-type: none"> ◆ present ◆ does not apply 	NA	present means the analyzed range contains off-scale dye sample peaks.

Searching the BioLIMS Database

Follow these steps to use the Collection Browser window to search the BioLIMS database for specific collections and fragments.

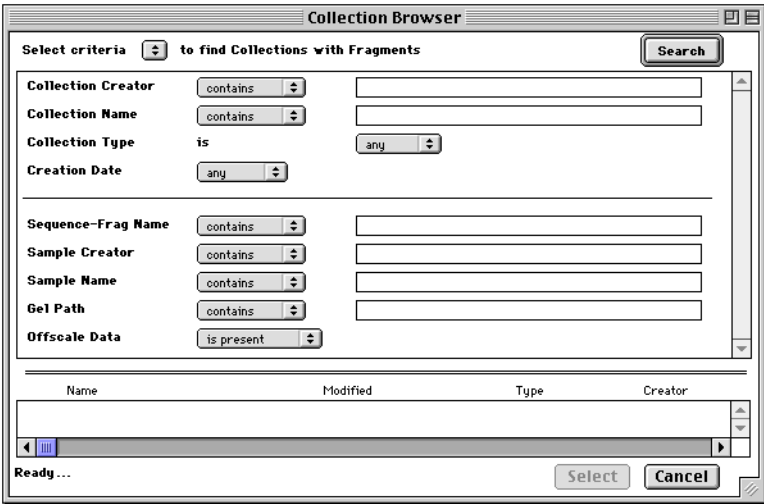
To search the BioLIMS database:

Step	Action
1	<p>Note From the Select Criteria pop-up menu, select the criteria by which you want to search</p> <p>Note To list all of the items in the BioLIMS database, perform the search with no criteria specified. For large databases, this process may be slow.</p>

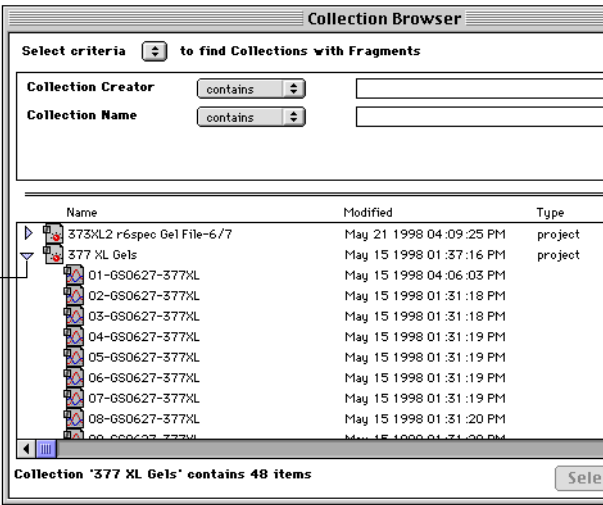


2	To use the pop-up menu:										
	<table border="1"> <thead> <tr> <th>Choose menu items...</th> <th>To define the search by...</th> <th>See page</th> </tr> </thead> <tbody> <tr> <td>above the horizontal line</td> <td>Collection Search Criteria</td> <td>E-24</td> </tr> <tr> <td>below the horizontal line</td> <td>Fragment Search Criteria</td> <td>E-25</td> </tr> </tbody> </table>	Choose menu items...	To define the search by...	See page	above the horizontal line	Collection Search Criteria	E-24	below the horizontal line	Fragment Search Criteria	E-25	
Choose menu items...	To define the search by...	See page									
above the horizontal line	Collection Search Criteria	E-24									
below the horizontal line	Fragment Search Criteria	E-25									
	<p>Note As you choose items from the pop-up menu, a black dot appears next to the item on the menu and the item is added to either the collection search criteria or the fragment name search criteria section of the window.</p>										

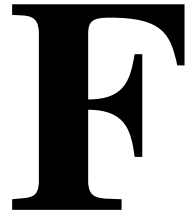
To search the BioLIMS database: *(continued)*

Step	Action
	<p>The following is an example of the Collection Browser window showing four collection search criteria and five fragment search criteria.</p>  <p>3 Use the pop-up menus and text fields to define your search query. Refer to “Collection Search Criteria” on page E-24, and “Fragment Search Criteria” on page E-25, for details about the search criteria. When you are satisfied with the search, click Search. The results of the search appear in the lower portion of the window.</p> <p>Note Collections returned by the Collection Browser window must match all of the collection criteria and contain at least one fragment that matches all of the fragment criteria.</p>

To search the BioLIMS database: *(continued)*

Step	Action
4	<p>To view the fragments contained in the collections, click the small triangle to the left of the collection name.</p>  <p>Small triangle</p>
5	Select fragments.

Troubleshooting the BioLIMS Database



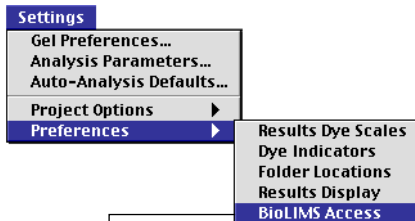
Introduction

In This Appendix Topics in this appendix include the following:

Topic	See Page
If the BioLIMS™ Preference Page Does Not Appear	F-2
About Troubleshooting the Client to Sybase Connection	F-3
Troubleshooting Process	F-15
Procedures for Troubleshooting the Client to Sybase Connection	F-6
About Troubleshooting the Client to Oracle Connection	F-14
Troubleshooting Process	F-15
Procedures for Troubleshooting the Client to Oracle Connection	F-17

If the BioLIMS™ Preference Page Does Not Appear

Problem If the BioLIMS™ Access page is not present after “Results Display” in the Preferences submenu, the GeneScan® Analysis Software is unable to find all the database support files and system extensions required to access the BioLIMS database.



This BioLIMS Access menu item will not appear if certain database support files are missing

Solution Be sure that all the Oracle® or Sybase SQL Server™ database support files are installed correctly.

If files are missing, reinstall the BioLIMS Client or Instrument package from the original CD-ROM disc.

About Troubleshooting the Client to Sybase Connection

Introduction A common source of difficulty using the BioLIMS System is establishing connection between the BioLIMS programs running on Macintosh® client computers and the BioLIMS database on the Sybase® Server.

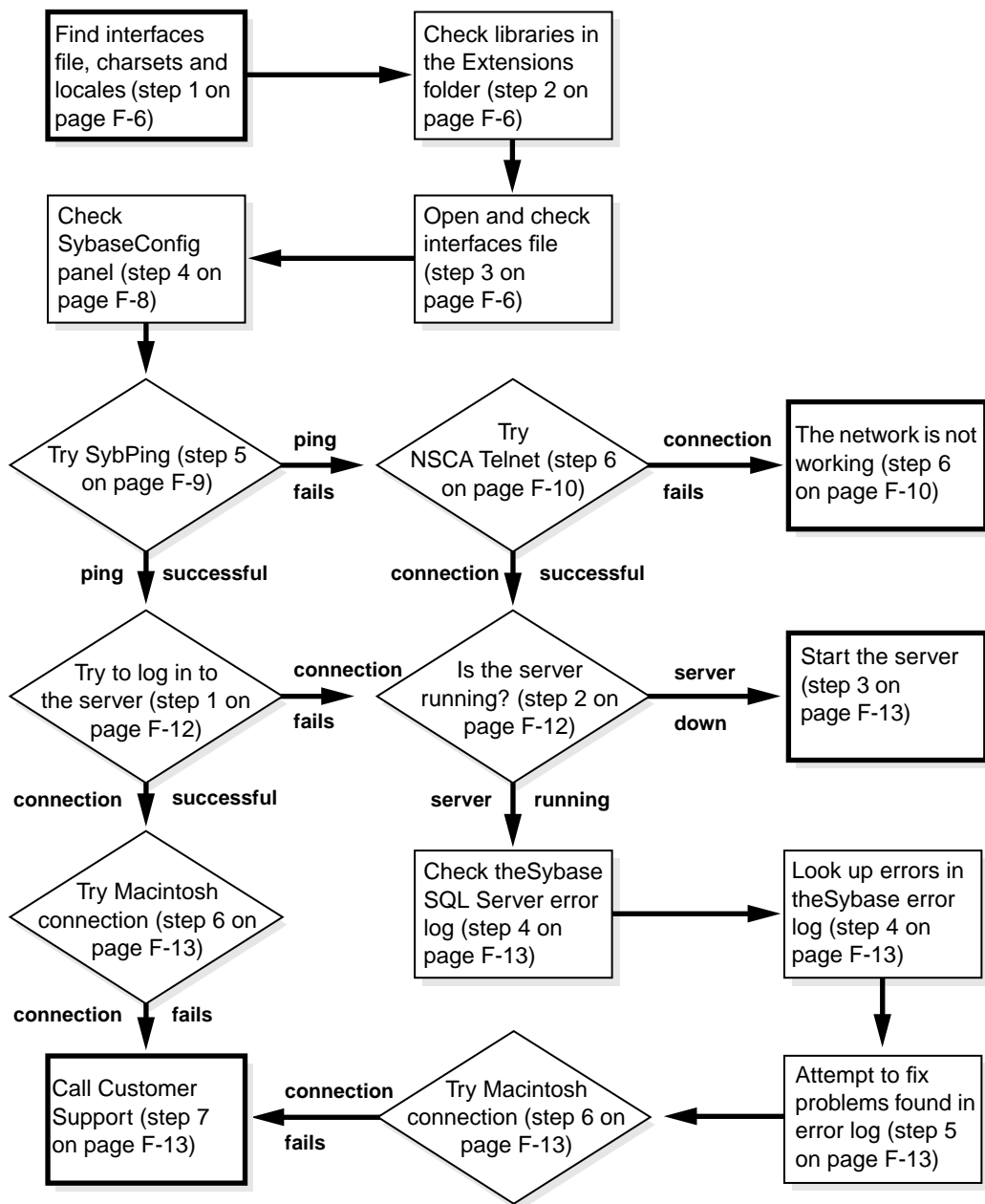
Note For information about troubleshooting an Oracle Server connection, see "About Troubleshooting the Client to Oracle Connection" on page F-14.

SybPing and Telnet Two programs have been provided to help with troubleshooting.

Program	Description
SybPing	A Sybase tool for testing network connections. (Look for this application in the BioLIMS:BioLIMS Extras:Sybase:bin folder.)
Telnet	<p>NCSA Telnet 2.6 is a program used for interactive access from a Macintosh client to a telnet host on TCP/IP networks. NCSA Telnet was developed by the National Center for Supercomputing Applications at the University of Illinois in Urbana/Champaign.</p> <p>NCSA Telnet 2.6 has been provided in the BioLIMS Extras folder.</p> <p>If you need help, their web site address is: http://www.ncsa.uiuc.edu/SDG/Software/Brochure/Overview/MacTelnet.overview.html</p> <p>You can also reach them at the address shown below. PE Applied Biosystems does not support NCSA Telnet.</p> <div style="border: 1px solid black; padding: 10px; margin: 10px 0;"> <p style="text-align: center; font-size: 2em; color: purple; font-family: cursive;">NCSA Telnet</p> <hr style="border: 0.5px solid black;"/> <p style="text-align: center; color: red;">Jim Browne Scott Bulmahn Jim Logan</p> <hr style="border: 0.5px solid black;"/> <p><i>Former Developers:</i> Tim Krauskopf, Geige Paulsen, Aaron Contorer, & Dave Whittington</p> <p><i>Contributors:</i> Bernt Budde, Basil Duval, Pascal Maes, Mark Tamsky, & Rick Watson</p> <p><i>Special Thanks:</i> The beta testers, the regulars on #macintosh and c.s.m.p., our tireless technical support team, and those who report the bugs.</p> <hr style="border: 0.5px solid black;"/> <p>NCSA Telnet is in the public domain. Comments and suggestions (catalog available) may be directed to:</p> <div style="display: flex; justify-content: space-between;"> <div style="width: 45%;"> <p>2.6 Think C 6.0 (March 1994)</p> </div> <div style="width: 45%; text-align: right;"> <p>NCSA Software Development 152 Computing Applications Bldg 605 E. Springfield Ave. Champaign, Ill 61820 mactelnet@ncsa.uiuc.edu</p> </div> </div> </div>

Troubleshooting Process

**Troubleshooting
Flow Chart** On the following page is a flow chart illustrating the process for troubleshooting database connection between the Macintosh Client and the Sybase Server. The step numbers given in the flow chart refer to the tables on pages F-6 to F-13 where the troubleshooting procedures are described in detail.



Procedures for Troubleshooting the Client to Sybase Connection

Introduction The procedure for troubleshooting the Client-Server connection is divided into two parts:

- ◆ Troubleshooting from the Macintosh Client (below).
- ◆ Troubleshooting from the Unix (Sybase) Server (page F-12).

Troubleshooting from the Macintosh Client

The step numbers in the following procedure correspond to the steps marked in the flow chart above.

IMPORTANT The documentation for each BioLIMS Macintosh computer program includes a section about setting up the database connection. Make sure you have followed that procedure carefully.

Troubleshooting the connection from the client (Macintosh computer):

Step	Action
1	Locate the interfaces file. Confirm that the folders charsets and locales are located in the same folder as the interfaces file. The default installation places interfaces, charsets, and locales in the BioLIMS:BioLIMS Extras:Sybase folder.
2	Confirm that the Sybase library files are installed into the Extensions folder in the System Folder. For more information, see "Required Sybase Extension Files" on page F-11. In particular, if the libtcp extension is missing or not turned on, the Sample 2DB Software crashes when connection to the database is attempted.
3	Open the interfaces file and confirm the server information is correct. An example of the Sybase-based BioLIMS SQL Server entry is shown below. <i>Sybase</i> query MacTCP mac_ether neuron.apldbio.com 2500 Note Refer to the table immediately below for an explanation of the BioLIMS Server entry.

Troubleshooting the connection from the client
(Macintosh computer): *(continued)*

Step	Action
The BioLIMS Server entry:	
Where...	Represents...
<i>SYBASE</i>	<p>the name you chose to call the BioLIMS Server. The server can have any name.</p> <p>For more information, see "About Server Names" on page E-17.</p>
query MacTCP mac_ether	<p>the part of the entry that is always the same.</p> <p>Note The tab preceding this phrase is required.</p>
neuron.apldbio. com	<p>the host and domain name of the Sybase-based BioLIMS SQL Server machine.</p> <p>In this example, neuron is the host name and apldbio.com is the domain name.</p> <p>You can also use an IP address. This information is available from your system administrator.</p>
2500	<p>the port number that the Sybase-based BioLIMS SQL Server is using to connect with the clients.</p> <p>This number is assigned to the server when it is installed.</p> <p>You can find the port number in the interfaces file, which is located in the home directory of the Sybase-based BioLIMS SQL Server.</p> <p>This is an example of the server entry in the UNIX server interfaces file:</p> <pre data-bbox="665 1104 1182 1218">## SYBASE on neuron ## Services: ## query tcp (2500) ...</pre> <p>The third line of this entry shows the port number (2500) of the BioLIMS Server.</p>

Troubleshooting the connection from the client
(Macintosh computer): *(continued)*

Step	Action												
4	Confirm that the SybaseConfig control panel is set up correctly:												
	<table border="1"> <thead> <tr> <th data-bbox="463 285 561 326">Step</th> <th data-bbox="561 285 1186 326">Action</th> </tr> </thead> <tbody> <tr> <td data-bbox="463 326 561 542">a.</td> <td data-bbox="561 326 1186 542"> <p>Open the SybaseConfig control panel.</p> <p>This control panel must be located in the Control Panels folder in the System folder.</p> <p>(If it is missing or disabled, the GeneScan Analysis Software crashes when connection to the database is attempted.)</p> </td> </tr> <tr> <td data-bbox="463 542 561 583">b.</td> <td data-bbox="561 542 1186 583">Click the Interfaces File button to open a browser box.</td> </tr> <tr> <td data-bbox="463 583 561 651">c.</td> <td data-bbox="561 583 1186 651">Use the browser box to locate and select the interfaces file that you located in step 1 above.</td> </tr> <tr> <td data-bbox="463 651 561 857">d.</td> <td data-bbox="561 651 1186 857"> <p>Confirm that the correct default server is chosen. Use the Default Server pop-up menu to select the default server.</p> <p>Note This is the same server you named for SYBASE in the BioLIMS Server entry table immediately above.</p> </td> </tr> <tr> <td data-bbox="463 857 561 928">e.</td> <td data-bbox="561 857 1186 928">Confirm that the Default Language pop-up menu is set to us_english.</td> </tr> </tbody> </table>	Step	Action	a.	<p>Open the SybaseConfig control panel.</p> <p>This control panel must be located in the Control Panels folder in the System folder.</p> <p>(If it is missing or disabled, the GeneScan Analysis Software crashes when connection to the database is attempted.)</p>	b.	Click the Interfaces File button to open a browser box.	c.	Use the browser box to locate and select the interfaces file that you located in step 1 above.	d.	<p>Confirm that the correct default server is chosen. Use the Default Server pop-up menu to select the default server.</p> <p>Note This is the same server you named for SYBASE in the BioLIMS Server entry table immediately above.</p>	e.	Confirm that the Default Language pop-up menu is set to us_english.
	Step	Action											
	a.	<p>Open the SybaseConfig control panel.</p> <p>This control panel must be located in the Control Panels folder in the System folder.</p> <p>(If it is missing or disabled, the GeneScan Analysis Software crashes when connection to the database is attempted.)</p>											
	b.	Click the Interfaces File button to open a browser box.											
	c.	Use the browser box to locate and select the interfaces file that you located in step 1 above.											
d.	<p>Confirm that the correct default server is chosen. Use the Default Server pop-up menu to select the default server.</p> <p>Note This is the same server you named for SYBASE in the BioLIMS Server entry table immediately above.</p>												
e.	Confirm that the Default Language pop-up menu is set to us_english.												

Troubleshooting the connection from the client
(Macintosh computer): *(continued)*

Step	Action														
5	<p data-bbox="517 215 1166 272">Use the program SybPing to confirm communication with the Sybase-based BioLIMS SQL Server.</p> <p data-bbox="517 289 1220 345">The default installation places SybPing into the BioLIMS: BioLIMS Extras: Sybase: bin folder.</p> <table border="1" data-bbox="517 391 1237 695" style="width: 100%; margin-top: 10px;"> <thead> <tr> <th data-bbox="517 391 611 425">Step</th> <th data-bbox="611 391 1237 425">Action</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 425 611 467" style="text-align: center;">a.</td> <td data-bbox="611 425 1237 467">Start the SybPing program.</td> </tr> <tr> <td data-bbox="517 467 611 581" style="text-align: center;">b.</td> <td data-bbox="611 467 1237 581"> Select the server from the Servers pop-up menu. (The servers shown in the pop-up menu are the servers listed in the interfaces file.) </td> </tr> <tr> <td data-bbox="517 581 611 695" style="text-align: center;">c.</td> <td data-bbox="611 581 1237 695"> Click Ping. The program responds with a message of whether or not the Ping was successful. </td> </tr> </tbody> </table> <table border="1" data-bbox="517 748 1237 898" style="width: 100%; margin-top: 10px;"> <thead> <tr> <th data-bbox="517 748 727 782">If the Ping is...</th> <th data-bbox="727 748 1237 782">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 782 727 857" style="text-align: center;">successful</td> <td data-bbox="727 782 1237 857">skip to "Troubleshooting the client connection from the Sybase SQL Server:" on page F-12.</td> </tr> <tr> <td data-bbox="517 857 727 898" style="text-align: center;">unsuccessful</td> <td data-bbox="727 857 1237 898">continue to step 6.</td> </tr> </tbody> </table>	Step	Action	a.	Start the SybPing program.	b.	Select the server from the Servers pop-up menu. (The servers shown in the pop-up menu are the servers listed in the interfaces file.)	c.	Click Ping. The program responds with a message of whether or not the Ping was successful.	If the Ping is...	Then...	successful	skip to "Troubleshooting the client connection from the Sybase SQL Server:" on page F-12.	unsuccessful	continue to step 6.
Step	Action														
a.	Start the SybPing program.														
b.	Select the server from the Servers pop-up menu. (The servers shown in the pop-up menu are the servers listed in the interfaces file.)														
c.	Click Ping. The program responds with a message of whether or not the Ping was successful.														
If the Ping is...	Then...														
successful	skip to "Troubleshooting the client connection from the Sybase SQL Server:" on page F-12.														
unsuccessful	continue to step 6.														

Troubleshooting the connection from the client
(Macintosh computer): *(continued)*

Step	Action								
6	If the Ping is unsuccessful: Open the program NCSA Telnet. The default installation places NCSA Telnet into the BioLIMS: BioLIMS Extras folder.								
	<table border="1"> <thead> <tr> <th>Step</th> <th>Action</th> </tr> </thead> <tbody> <tr> <td>a.</td> <td>Select Open Connection... from the File menu.</td> </tr> <tr> <td>b.</td> <td>Enter the host name (from the interfaces file) into the Host/Session Name text field.</td> </tr> <tr> <td>c.</td> <td>Click Connect.</td> </tr> </tbody> </table>	Step	Action	a.	Select Open Connection... from the File menu.	b.	Enter the host name (from the interfaces file) into the Host/Session Name text field.	c.	Click Connect.
	Step	Action							
	a.	Select Open Connection... from the File menu.							
	b.	Enter the host name (from the interfaces file) into the Host/Session Name text field.							
	c.	Click Connect.							
	<table border="1"> <thead> <tr> <th>If the connection is...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>successful</td> <td>a window is displayed with a UNIX login prompt. Skip to step 2 of "Troubleshooting the client connection from the Sybase SQL Server:" on page F-12.</td> </tr> <tr> <td>unsuccessful</td> <td>an error message is displayed. The network is not working. Call your network administrator. Note If the host.domain name does not work, try the IP address of the host. This number can be found in the /etc/hosts file on the UNIX Server.</td> </tr> </tbody> </table>	If the connection is...	Then...	successful	a window is displayed with a UNIX login prompt. Skip to step 2 of "Troubleshooting the client connection from the Sybase SQL Server:" on page F-12.	unsuccessful	an error message is displayed. The network is not working. Call your network administrator. Note If the host.domain name does not work, try the IP address of the host. This number can be found in the /etc/hosts file on the UNIX Server.		
	If the connection is...	Then...							
	successful	a window is displayed with a UNIX login prompt. Skip to step 2 of "Troubleshooting the client connection from the Sybase SQL Server:" on page F-12.							
	unsuccessful	an error message is displayed. The network is not working. Call your network administrator. Note If the host.domain name does not work, try the IP address of the host. This number can be found in the /etc/hosts file on the UNIX Server.							

**Required Sybase
Extension Files**

The following Sybase library files are placed in the Extensions folder in the System folder by the BioLIMS Client and instrument installers. These files are required for connection to the BioLIMS database.

libblk	libintl
libcomn	libsybdb
libcs	libtcl
libct	libtcp
libctb	

**Required Oracle
Extension Files**

The following Oracle library files are placed in the Extensions folder in the System folder by the BioLIMS Client and instrument installers. These files are required for connection to the BioLIMS database.

OracleCore23Lib	OracleOra71Lib
OracleKernel71Lib	OraclePlsql21Lib
OracleNetNLLib	OraclePstd21Lib
OracleNetTCPLib	OracleRuntime13Lib
OracleNetTNSLib	OracleSql16Lib
OracleNetTNSTCPLib	OracleTNSATKLib
OracleNLS23Lib	OracleVsoci71Lib
OracleOci71Lib	

Troubleshooting from the Unix (Sybase) Server

The step numbers in the following procedure correspond to the steps marked in the flow chart on page F-5. Step 1 below corresponds to the diamond "Try to log in to the server" on page F-5.

You need to have the account name and password for the Sybase user on the UNIX system that runs the Sybase SQL Server. If you do not have access to the Sybase user account, you should ask your database administrator to carry out the following procedure

Troubleshooting the client connection from the Sybase SQL Server:

Step	Action										
1	Log in to the Sybase user account on the UNIX Server.										
	Try to connect to the Sybase SQL Server with isql.										
	Use the same client user name as the one entered in the BioLIMS access dialog box.										
	For example:										
	<pre>% isql -U george -P george1. -S SYBASE 1> use sfdb 2> go 1> quit</pre>										
	<table border="1"> <thead> <tr> <th>Where...</th> <th>Represents the...</th> </tr> </thead> <tbody> <tr> <td>george</td> <td>client user name.</td> </tr> <tr> <td>george1</td> <td>client user password.</td> </tr> <tr> <td>SYBASE</td> <td>server name.</td> </tr> <tr> <td>sfdb</td> <td>name of the BioLIMS database.</td> </tr> </tbody> </table>	Where...	Represents the...	george	client user name.	george1	client user password.	SYBASE	server name.	sfdb	name of the BioLIMS database.
Where...	Represents the...										
george	client user name.										
george1	client user password.										
SYBASE	server name.										
sfdb	name of the BioLIMS database.										
	<table border="1"> <thead> <tr> <th>If the Login is...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>successful</td> <td>skip to step 6.</td> </tr> <tr> <td>unsuccessful</td> <td>continue to step 2.</td> </tr> </tbody> </table>	If the Login is...	Then...	successful	skip to step 6.	unsuccessful	continue to step 2.				
If the Login is...	Then...										
successful	skip to step 6.										
unsuccessful	continue to step 2.										
2	Use the show server script to find out if the Sybase SQL Server is running.										
	<table border="1"> <thead> <tr> <th>If the server is...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>running</td> <td>skip to step 4.</td> </tr> <tr> <td>not running</td> <td>continue to step 3.</td> </tr> </tbody> </table>	If the server is...	Then...	running	skip to step 4.	not running	continue to step 3.				
	If the server is...	Then...									
running	skip to step 4.										
not running	continue to step 3.										

Troubleshooting the client connection from the **Sybase SQL**

Step	Action						
3	<p>Ask your database administrator to restart the Sybase SQL Server.</p> <p>After restarting the server, try to connect to the database using isql as in step 1 on page F-12.</p> <p>If the connection still fails, go to step 7.</p>						
4	<p>Check the Sybase error log and make a note of any error messages.</p> <p>(The error log can be found in the install directory of the Sybase home directory and error messages are preceded by the string Msg.)</p>						
5	<p>Refer to the Sybooks documentation for an explanation of the SQL Server Error Messages. Attempt to fix the problem following the Sybook instructions.</p> <p>Try to connect to the database using isql as in step 1 on page F-12.</p> <table border="1" data-bbox="516 685 1233 805"> <thead> <tr> <th>If the connection is...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>successful</td> <td>continue to step 6.</td> </tr> <tr> <td>unsuccessful</td> <td>skip to step 7.</td> </tr> </tbody> </table>	If the connection is...	Then...	successful	continue to step 6.	unsuccessful	skip to step 7.
If the connection is...	Then...						
successful	continue to step 6.						
unsuccessful	skip to step 7.						
6	<p>Try once more to connect to the BioLIMS database from the BioLIMS Macintosh program.</p> <p>Make sure that the user name, password, database and server names are all typed correctly and are in the correct case.</p> <p>If the connection still fails, go to step 7.</p>						
7	<p>Call Customer Support. See "Technical Support" on page 1-20.</p>						

About Troubleshooting the Client to Oracle Connection

Introduction A common source of difficulty using the BioLIMS System is establishing connection between the BioLIMS programs running on Macintosh client computers and the BioLIMS database on an Oracle Server.

Note For information about troubleshooting a Sybase SQL Server connection, see "About Troubleshooting the Client to Sybase Connection" on page F-3.

Telnet To help with troubleshooting, NCSA Telnet 2.6 has been provided in the BioLIMS Extras folder.

NCSA Telnet 2.6 is a program used for interactive access from a Macintosh client to a telnet host on TCP/IP networks. NCSA Telnet was developed by the National Center for Supercomputing Applications at the University of Illinois in Urbana/Champaign.

If you need help, their web site address is:

<http://www.ncsa.uiuc.edu/SDG/Software/Brochure/Overview/MacTelnet.overview.html>

You can also reach them at the address shown below.

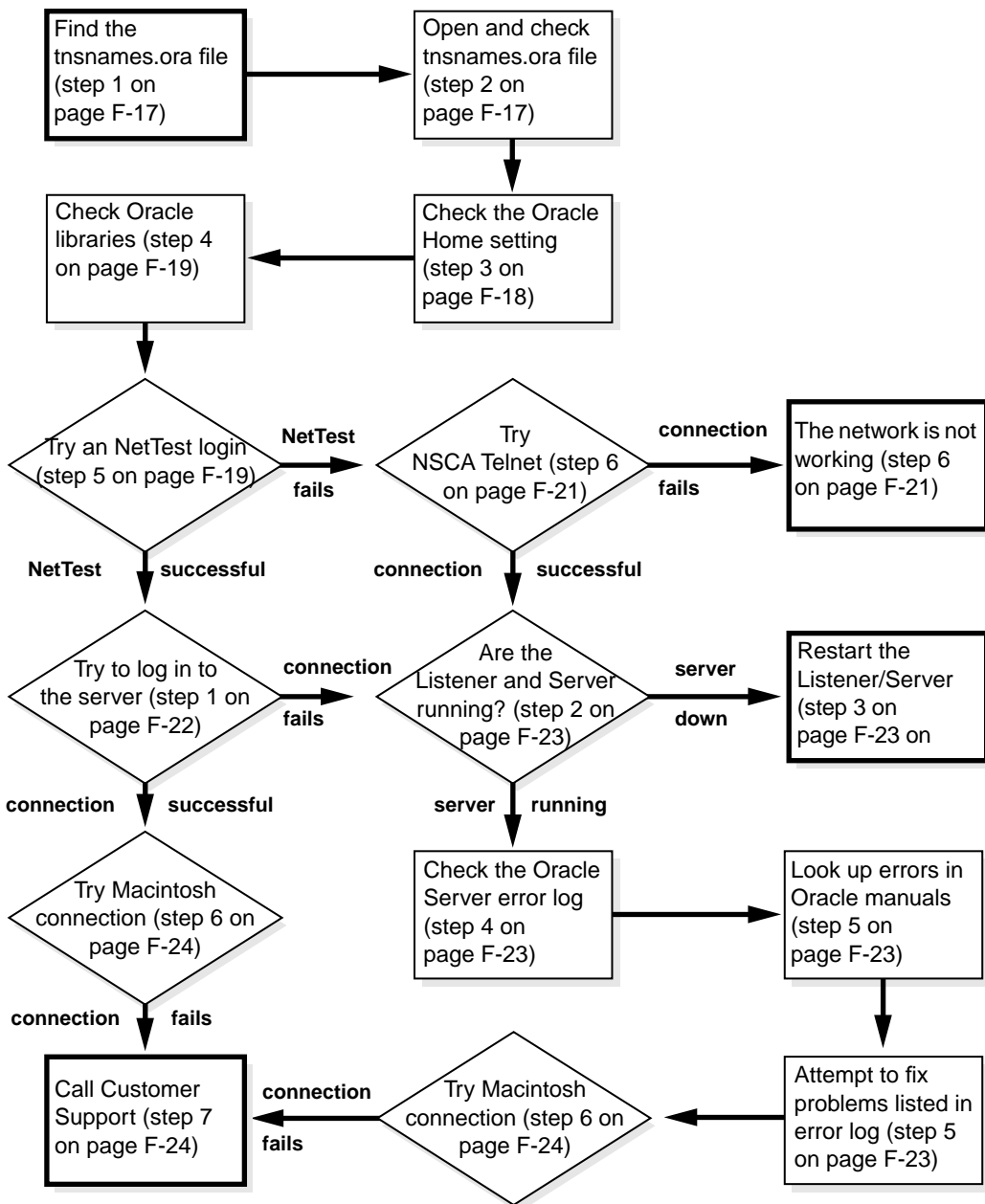
PE Applied Biosystems does not support NCSA Telnet.

NCSA Telnet		
Jim Browne	Scott Bulmahn	Jim Logan
<i>Former Developers:</i> Tim Krauskopf, Gaige Paulsen, Aaron Contorer, & Dave Whittington		
<i>Contributors:</i> Bernt Budde, Basil Duval, Pascal Maes, Mark Tamsky, & Rick Watson		
<i>Special Thanks:</i> The beta testers, the regulars on #macintosh and c.s.m.p., our tireless technical support team, and those who report the bugs.		
NCSA Telnet is in the public domain. Comments and suggestions (catalog available) may be directed to:		NCSA Software Development 152 Computing Applications Bldg 605 E. Springfield Ave. Champaign, Ill 61820 mactelnet@ncsa.uiuc.edu
2.6 Think C 6.0 (March 1994)		

Troubleshooting Process

Troubleshooting Flow Chart

On the following page is a flow chart illustrating the process for troubleshooting the database connection between the Macintosh Client and the Oracle Server. The step numbers given in the flow chart refer to the tables on pages F-17 to F-22 where the troubleshooting procedures are described in detail.



Procedures for Troubleshooting the Client to Oracle Connection

Introduction The procedure for troubleshooting the Client-Server connection is divided into two parts:

- ◆ Troubleshooting from the Macintosh Client (below).
- ◆ Troubleshooting from the Unix (Oracle) Server (page F-22).

Troubleshooting from the Macintosh Client The step numbers in the following procedure correspond to the steps marked in the flow chart above.

IMPORTANT The documentation for each BioLIMS Macintosh program includes a section about setting up the database connection. Ensure that you have followed that procedure carefully.

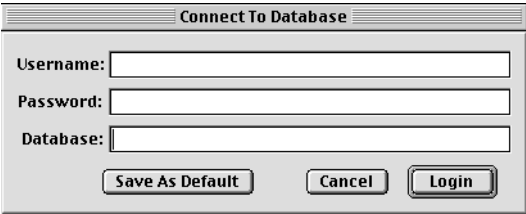
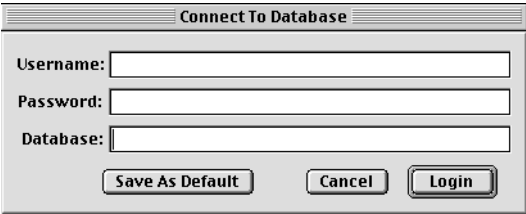
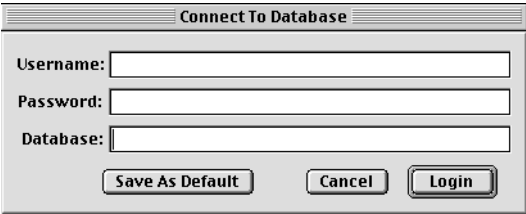
Troubleshooting the connection from the client (Macintosh)

Step	Action
1	Locate the tnsnames.ora file. The default installation places the tnsnames.ora file in the BioLIMS 2.0:BioLIMS Extras:Oracle:Network:Admin folder.
2	Open the tnsnames.ora file and confirm that the server information is correct. An example of the Oracle-based BioLIMS Server entry is shown below.
<pre>Oramozart = (DESCRIPTION= (ADDRESS= (PROTOCOL=TCP) (host=mozart) (port=1521)) (CONNECT_DATA=(SID=WG733)))</pre>	
Note	Refer to the table immediately below for an explanation of the BioLIMS Server entry.

Troubleshooting the connection from the client (Macintosh) *(continued)*

Step	Action
The BioLIMS Server entry:	
Where...	Represents...
<i>Oramozart</i>	<p>the name you chose to call the Oracle-based BioLIMS database connection.</p> <p>This can have any name.</p> <p>See "About Server Names" on page E-17</p>
<i>mozart</i>	<p>the host of the Oracle-based BioLIMS Server.</p> <p>In this example, mozart is the host name.</p> <p>You can also use a host and domain name such as mozart.apldbio.com or an IP address such as 167.116.100.61.</p> <p>This information is available from your system administrator. The IP address can also be found in the /etc/hosts file on the UNIX Server.</p>
1521	<p>the port number that the Oracle-based BioLIMS Server is using to connect with the clients.</p> <p>This number is assigned to the server when it is installed.</p> <p>You can find the port number in the tnsnames.ora file that is located in the \$ORACLE_HOME/network/admin directory on the Oracle-based BioLIMS Server.</p>
WG733	<p>the Oracle database system identifier (sid) defined at server installation.</p> <p>You can find the port number in the tnsnames.ora file that is located in the \$ORACLE_HOME/network/admin directory on the Oracle-based BioLIMS Server.</p>
3	<p>Confirm that the Oracle Home is set correctly:</p> <ol style="list-style-type: none"> a. Open the Set Oracle Home program. At installation this program is placed into the BioLIMS 2.0:BioLIMS Extras:Oracle:Applications folder. b. Use the file browser to locate and select the Oracle folder.

Troubleshooting the connection from the client (Macintosh) *(continued)*

Step	Action												
4	<p>Confirm that the Oracle library files are installed.</p> <p>See “Required Oracle Extension Files” on page F-11.</p>												
5	<p>Use the program NetTest to confirm communication with the Oracle Server.</p> <p>At installation, NetTest is placed into the BioLIMS 2.0:BioLIMS Extras:Oracle:Applications:Networking folder.</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th>Step</th> <th>Action</th> </tr> </thead> <tbody> <tr> <td>a.</td> <td> <p>Open the NetTest program.</p> <p>Note No window appears for the NetTest program, but the menu bar changes.</p> </td> </tr> <tr> <td>b.</td> <td> <p>From the Database menu, choose Logon.</p> <p>The Connect To Database dialog box appears.</p> <div style="text-align: center;">  </div> </td> </tr> <tr> <td>c.</td> <td> <p>d. Enter the Username and Password for the database.</p> <p>Note These are the same Username and Password you use to log into BioLIMS (e.g., george/george1).</p> </td> </tr> <tr> <td>e.</td> <td> <p>In the Database field, enter the alias name of the Oracle Server from the tnsnames.ora file (e.g., Oramozart).</p> </td> </tr> <tr> <td>f.</td> <td> <p>Click Login.</p> <p>The message “Attempting Connection...” appears, then the Result Explanation dialog box appears.</p> </td> </tr> </tbody> </table>	Step	Action	a.	<p>Open the NetTest program.</p> <p>Note No window appears for the NetTest program, but the menu bar changes.</p>	b.	<p>From the Database menu, choose Logon.</p> <p>The Connect To Database dialog box appears.</p> <div style="text-align: center;">  </div>	c.	<p>d. Enter the Username and Password for the database.</p> <p>Note These are the same Username and Password you use to log into BioLIMS (e.g., george/george1).</p>	e.	<p>In the Database field, enter the alias name of the Oracle Server from the tnsnames.ora file (e.g., Oramozart).</p>	f.	<p>Click Login.</p> <p>The message “Attempting Connection...” appears, then the Result Explanation dialog box appears.</p>
Step	Action												
a.	<p>Open the NetTest program.</p> <p>Note No window appears for the NetTest program, but the menu bar changes.</p>												
b.	<p>From the Database menu, choose Logon.</p> <p>The Connect To Database dialog box appears.</p> <div style="text-align: center;">  </div>												
c.	<p>d. Enter the Username and Password for the database.</p> <p>Note These are the same Username and Password you use to log into BioLIMS (e.g., george/george1).</p>												
e.	<p>In the Database field, enter the alias name of the Oracle Server from the tnsnames.ora file (e.g., Oramozart).</p>												
f.	<p>Click Login.</p> <p>The message “Attempting Connection...” appears, then the Result Explanation dialog box appears.</p>												

Troubleshooting the connection from the client (Macintosh) *(continued)*

Step	Action						
	<table border="1"> <thead> <tr> <th data-bbox="471 207 749 246">If the login is...</th> <th data-bbox="749 207 1176 246">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="471 246 749 634">successful</td> <td data-bbox="749 246 1176 634"> <p>the Result Explanation dialog box displays the following:</p> <ul style="list-style-type: none"> ◆ Result field: "ORA-00000: normal, successful completion." ◆ Connected To field: Oracle Server version information. ◆ Comments field: "Your SQL*Net setup appears to be in working order." <p>Skip to "Troubleshooting from the Unix (Oracle) Server" on page F-22.</p> </td> </tr> <tr> <td data-bbox="471 634 749 889">unsuccessful</td> <td data-bbox="749 634 1176 889"> <p>the Result Explanation dialog box displays the following:</p> <ul style="list-style-type: none"> ◆ Result field: the Oracle error encountered. ◆ Comments field: A detailed explanation of the login failure. <p>Continue to step 6.</p> </td> </tr> </tbody> </table>	If the login is...	Then...	successful	<p>the Result Explanation dialog box displays the following:</p> <ul style="list-style-type: none"> ◆ Result field: "ORA-00000: normal, successful completion." ◆ Connected To field: Oracle Server version information. ◆ Comments field: "Your SQL*Net setup appears to be in working order." <p>Skip to "Troubleshooting from the Unix (Oracle) Server" on page F-22.</p>	unsuccessful	<p>the Result Explanation dialog box displays the following:</p> <ul style="list-style-type: none"> ◆ Result field: the Oracle error encountered. ◆ Comments field: A detailed explanation of the login failure. <p>Continue to step 6.</p>
If the login is...	Then...						
successful	<p>the Result Explanation dialog box displays the following:</p> <ul style="list-style-type: none"> ◆ Result field: "ORA-00000: normal, successful completion." ◆ Connected To field: Oracle Server version information. ◆ Comments field: "Your SQL*Net setup appears to be in working order." <p>Skip to "Troubleshooting from the Unix (Oracle) Server" on page F-22.</p>						
unsuccessful	<p>the Result Explanation dialog box displays the following:</p> <ul style="list-style-type: none"> ◆ Result field: the Oracle error encountered. ◆ Comments field: A detailed explanation of the login failure. <p>Continue to step 6.</p>						

Troubleshooting the connection from the client (Macintosh) *(continued)*

Step	Action														
6	<p data-bbox="517 183 893 207">If the NetTest login is unsuccessful:</p> <p data-bbox="517 228 1229 280">Open the program NCSA Telnet. The default installer places NCSA Telnet into the BioLIMS 2.0:BioLIMS Extras folder.</p> <table border="1" data-bbox="517 329 1236 621" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th data-bbox="517 329 611 370">Step</th> <th data-bbox="611 329 1236 370">Action</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 370 611 410" style="text-align: center;">a.</td> <td data-bbox="611 370 1236 410">Select Open Connection... from the File menu.</td> </tr> <tr> <td data-bbox="517 410 611 581" style="text-align: center;">b.</td> <td data-bbox="611 410 1236 581"> Enter the host name (from the tnsnames.ora file) into the Host/Session Name text field. Note You can enter the host, host and domain name, or IP address here. See step 2, "mozart," on page F-18 for more information. </td> </tr> <tr> <td data-bbox="517 581 611 621" style="text-align: center;">c.</td> <td data-bbox="611 581 1236 621">Click Connect.</td> </tr> </tbody> </table> <table border="1" data-bbox="517 678 1236 1226" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th data-bbox="517 678 786 719">If the connection is...</th> <th data-bbox="786 678 1236 719">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="517 719 786 889" style="text-align: center;">successful</td> <td data-bbox="786 719 1236 889"> a window is displayed with a Unix login prompt. Skip to step 2 on page F-12 of "Troubleshooting from the Unix (Oracle) Server." </td> </tr> <tr> <td data-bbox="517 889 786 1226" style="text-align: center;">unsuccessful</td> <td data-bbox="786 889 1236 1226"> an error message is displayed. The network is not working. Call your network administrator. Note If you entered the host name in step b above and the connection fails, try entering the host.domain name or the IP address before calling your network administrator. See step 2, "mozart," on page F-18 for more information. </td> </tr> </tbody> </table>	Step	Action	a.	Select Open Connection... from the File menu.	b.	Enter the host name (from the tnsnames.ora file) into the Host/Session Name text field. Note You can enter the host, host and domain name, or IP address here. See step 2, "mozart," on page F-18 for more information.	c.	Click Connect.	If the connection is...	Then...	successful	a window is displayed with a Unix login prompt. Skip to step 2 on page F-12 of "Troubleshooting from the Unix (Oracle) Server."	unsuccessful	an error message is displayed. The network is not working. Call your network administrator. Note If you entered the host name in step b above and the connection fails, try entering the host.domain name or the IP address before calling your network administrator. See step 2, "mozart," on page F-18 for more information.
Step	Action														
a.	Select Open Connection... from the File menu.														
b.	Enter the host name (from the tnsnames.ora file) into the Host/Session Name text field. Note You can enter the host, host and domain name, or IP address here. See step 2, "mozart," on page F-18 for more information.														
c.	Click Connect.														
If the connection is...	Then...														
successful	a window is displayed with a Unix login prompt. Skip to step 2 on page F-12 of "Troubleshooting from the Unix (Oracle) Server."														
unsuccessful	an error message is displayed. The network is not working. Call your network administrator. Note If you entered the host name in step b above and the connection fails, try entering the host.domain name or the IP address before calling your network administrator. See step 2, "mozart," on page F-18 for more information.														

Troubleshooting from the Unix (Oracle) Server

The step numbers in the following procedure correspond to the steps marked in the flow chart on page F-16.

You need to have the account name and password for the Oracle user on the UNIX system that runs the Oracle Server. If you do not have access to the Oracle user account, you should ask your database administrator to carry out the following procedure.

Troubleshooting the client connection from the Oracle Server:

Step	Action												
1	<p>Log in to the Oracle user account on the UNIX Server.</p> <p>Try to connect to the Oracle Server with SQL*Plus.</p> <p>Use the same client user name and password as the one entered in the BioLIMS access dialog box. For example:</p> <pre>% sqlplus george/george1 SQL>exit</pre> <table border="1" data-bbox="470 719 1189 837"> <thead> <tr> <th>Where...</th> <th>Represents the...</th> </tr> </thead> <tbody> <tr> <td>george</td> <td>client user name.</td> </tr> <tr> <td>george1</td> <td>client user password.</td> </tr> </tbody> </table> <table border="1" data-bbox="470 889 1189 1008"> <thead> <tr> <th>If the Login is...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>successful</td> <td>skip to step 6.</td> </tr> <tr> <td>unsuccessful</td> <td>continue to step 2.</td> </tr> </tbody> </table>	Where...	Represents the...	george	client user name.	george1	client user password.	If the Login is...	Then...	successful	skip to step 6.	unsuccessful	continue to step 2.
Where...	Represents the...												
george	client user name.												
george1	client user password.												
If the Login is...	Then...												
successful	skip to step 6.												
unsuccessful	continue to step 2.												

Troubleshooting the client connection from the Oracle Server: *(continued)*

Step	Action												
2	<p>Find out if the Listener Process is running.</p> <p>As the Oracle user, type:</p> <pre>% lsnrctl status</pre> <table border="1" data-bbox="514 354 1239 824"> <thead> <tr> <th data-bbox="514 354 705 423">If the Listener Process is...</th> <th data-bbox="705 354 1239 423">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="514 423 705 461">not running</td> <td data-bbox="705 423 1239 461">continue to step 3.</td> </tr> <tr> <td data-bbox="514 461 705 634">running</td> <td data-bbox="705 461 1239 634"> <p>check that the server is running.</p> <p>Type:</p> <pre>% ps -ef grep ora_</pre> <table border="1" data-bbox="713 634 1231 824"> <thead> <tr> <th data-bbox="713 634 974 704">If server processes are...</th> <th data-bbox="974 634 1231 704">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="713 704 974 774">displayed</td> <td data-bbox="974 704 1231 774">server is running. Skip to step 4.</td> </tr> <tr> <td data-bbox="713 774 974 824">not displayed</td> <td data-bbox="974 774 1231 824">continue to step 3.</td> </tr> </tbody> </table> </td> </tr> </tbody> </table>	If the Listener Process is...	Then...	not running	continue to step 3.	running	<p>check that the server is running.</p> <p>Type:</p> <pre>% ps -ef grep ora_</pre> <table border="1" data-bbox="713 634 1231 824"> <thead> <tr> <th data-bbox="713 634 974 704">If server processes are...</th> <th data-bbox="974 634 1231 704">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="713 704 974 774">displayed</td> <td data-bbox="974 704 1231 774">server is running. Skip to step 4.</td> </tr> <tr> <td data-bbox="713 774 974 824">not displayed</td> <td data-bbox="974 774 1231 824">continue to step 3.</td> </tr> </tbody> </table>	If server processes are...	Then...	displayed	server is running. Skip to step 4.	not displayed	continue to step 3.
If the Listener Process is...	Then...												
not running	continue to step 3.												
running	<p>check that the server is running.</p> <p>Type:</p> <pre>% ps -ef grep ora_</pre> <table border="1" data-bbox="713 634 1231 824"> <thead> <tr> <th data-bbox="713 634 974 704">If server processes are...</th> <th data-bbox="974 634 1231 704">Then...</th> </tr> </thead> <tbody> <tr> <td data-bbox="713 704 974 774">displayed</td> <td data-bbox="974 704 1231 774">server is running. Skip to step 4.</td> </tr> <tr> <td data-bbox="713 774 974 824">not displayed</td> <td data-bbox="974 774 1231 824">continue to step 3.</td> </tr> </tbody> </table>	If server processes are...	Then...	displayed	server is running. Skip to step 4.	not displayed	continue to step 3.						
If server processes are...	Then...												
displayed	server is running. Skip to step 4.												
not displayed	continue to step 3.												
3	<p>See your system administrator for help restarting the Listener Process or Server.</p> <p>After restarting, try to connect to the database using SQL*Plus as in step 1 on page F-22.</p> <p>If the connection still fails, go to step 7.</p>												
4	<p>Check the Oracle error log and make a note of any error messages.</p> <p>(The error log can be found in the install directory of the Oracle home directory and error messages are preceded by the string Msg.)</p> <p>Note Refer to the Oracle documentation for an explanation of the error messages.</p>												
5	<p>Attempt to fix the problem following the documentation instructions.</p> <p>Try to connect to the database using SQL*Plus as in step 1 on page F-22.</p> <p>If the connection still fails, go to step 7.</p>												

Troubleshooting the client connection from the Oracle Server: *(continued)*

Step	Action
6	Try once more to connect to the BioLIMS database from the BioLIMS Macintosh program. Ensure that the user name, password, database, and server names are all typed correctly and are in the correct case. If the connection still fails, go to step 7.
7	Call Customer Support. See "Technical Support" on page 1-20.

License and Warranty



Software License and Warranty

**Perkin-Elmer
Software License
and Limited
Product Warranty**

PURCHASER, CAREFULLY READ THE FOLLOWING TERMS AND CONDITIONS (THE "AGREEMENT"), WHICH APPLY TO THE SOFTWARE ENCLOSED (THE "SOFTWARE").

YOUR OPENING OF THIS PACKAGE INDICATES YOUR ACCEPTANCE OF THESE TERMS AND CONDITIONS. IF YOU DO NOT ACCEPT THEM, PROMPTLY RETURN THE COMPLETE PACKAGE AND YOUR MONEY WILL BE RETURNED. THE LAW PROVIDES FOR CIVIL AND CRIMINAL PENALTIES FOR ANYONE WHO VIOLATES THE LAWS OF COPYRIGHT.

Copyright

The SOFTWARE, including its structure, organization, code, user interface, and associated documentation, is a proprietary product of The Perkin-Elmer Corporation and is protected by international laws of copyright. Title to the SOFTWARE, and to any and all portion(s) of the SOFTWARE shall at all times remain with Perkin-Elmer.

License

1. You may use the SOFTWARE on a single computer (or on a single network, if your software is designated as a network version). You may transfer the SOFTWARE to another single computer (or network, if a network version), so long as you first delete the SOFTWARE from the previous computer or network. You may never have operational SOFTWARE on more than one computer (or more than one network, if a network version) per original copy of the SOFTWARE at any time.
2. You may make one copy of the SOFTWARE for backup purposes.
3. You may transfer the SOFTWARE to another party, but only if the other party agrees in writing with Perkin-Elmer to accept the terms and conditions of this Agreement. If you transfer the SOFTWARE to another

party, you must immediately transfer all copies to that party, or destroy those not transferred. Any such transfer terminates your license.

-
- Restrictions**
1. You may not copy, transfer, rent, modify, use, or merge the SOFTWARE, or the associated documentation, in whole or in part, except as expressly permitted in this Agreement.
 2. You may not reverse assemble, decompile, or otherwise reverse engineer the SOFTWARE.

Limited Warranty

For a period of 90 days after purchase of the SOFTWARE, Perkin-Elmer warrants that the SOFTWARE will function substantially as described in the documentation supplied by Perkin-Elmer with the SOFTWARE. If you discover an error which causes substantial deviation from that documentation, send a written notification to Perkin-Elmer. Upon receiving such notification, if Perkin-Elmer is able to reliably reproduce that error at its facility, then Perkin-Elmer will do one of the following at its sole option: (i) correct the error in a subsequent release of the SOFTWARE, which shall be supplied to you free of charge, or (ii) accept a return of the SOFTWARE from you, and refund the purchase price received for the SOFTWARE. Perkin-Elmer does not warrant that the SOFTWARE will meet your requirements, will be error-free, or will conform exactly to the documentation. Any sample or model used in connection with this Agreement is for illustrative purposes only, is not part of the basis of the bargain, and is not to be construed as a warranty that the SOFTWARE will conform to the sample or model.

Limitation Of Liability

EXCEPT AS SPECIFICALLY STATED IN THIS AGREEMENT, THE SOFTWARE IS PROVIDED AND LICENSED "AS IS". THE ABOVE WARRANTY IS GIVEN IN LIEU OF ALL OTHER WARRANTIES, EXPRESSED OR IMPLIED, INCLUDING THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. NOTWITHSTANDING ANY FAILURE OF THE CENTRAL PURPOSE OF ANY LIMITED REMEDY, PERKIN-ELMER'S LIABILITY FOR BREACH OF WARRANTY SHALL BE LIMITED TO A REFUND OF THE PURCHASE PRICE FOR SUCH PRODUCT. IN NO EVENT WILL PERKIN-ELMER BE LIABLE FOR ANY OTHER DAMAGES, INCLUDING INCIDENTAL OR CONSEQUENTIAL DAMAGES, EVEN IF PERKIN-ELMER HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

Term You may terminate this Agreement by destroying all copies of the SOFTWARE and documentation. Perkin-Elmer may terminate this Agreement if you fail to comply with any or all of its terms, in which case you agree to return to Perkin-Elmer all copies of the SOFTWARE and associated documentation.

-
-
- Miscellaneous**
1. Failure to enforce any of the terms and conditions of this Agreement by either party shall not be deemed a waiver of any rights and privileges under this Agreement.
 2. In case any one or more of the provisions of this Agreement for any reason shall be held to be invalid, illegal, or unenforceable in any respect, such invalidity, illegality, or unenforceability shall not affect any other provisions of this Agreement, and this Agreement shall be construed as if such invalid, illegal, or unenforceable provisions had never been contained herein.
 3. This Agreement shall be construed and governed by the laws of the State of California.
 4. This Agreement and the Perkin-Elmer Sales Quotation constitute the entire agreement between Perkin-Elmer and you concerning the SOFTWARE.
-
-

Glossary

This glossary defines special terminology used in the *GeneScan Analysis Software User's Manual*. The terms are listed in alphabetical order. Many terms are defined in the text of the manual, so if you do not find a term here, check the index to see if you can locate it in the manual.

align-by-size curve Curve created by the GeneScan® Analysis Software for aligning data by size. The software calculates a best-fit, least-squares curve for all samples. This is a third-order curve when you use the Third Order Least Squares size calling method; for all other size calling methods it is a second-order curve. This curve is black in the Standard Sizing Curve window, although when the sizing curve and this curve match, they overlap so you see only the sizing curve. See also “size calling curve,” *and* “size standard spline interpolation curve.”

analysis parameters Options that specify certain ranges and methods used during analysis using the GeneScan Analysis Software. The software has default analysis parameters that are stored in the project itself. These parameters apply globally, unless you create your own parameters files for use with specific protocols. See “Defining Analysis Parameters” on page 5-18.

baselining Adjusting the baselines of detected dye colors to the same level for a better comparison of relative signal intensity.

channels (ABI™ 373 and ABI PRISM® 377 instruments) Theoretical divisions across the read region of a gel where the Data Collection software samples the data. The number of wells in the loading comb used determines the approximate number of channels assigned per lane of the gel (for instance, with a 36-well comb, one lane is approximately five channels). When the GeneScan Analysis Software tracks a gel, it places the tracker line for each lane in the channel showing the strongest fluorescent signal. The data from that channel and the adjacent channels on each side are averaged to determine the raw data for the Sample file. You set the number of channels to be averaged in the Gel Processing Parameters.

data point The ABI 373 and ABI PRISM 377 Data Collection software samples data 194 times as it scans across the gel (194 or 388 times for the XL upgrades). The ABI PRISM® 310 Data Collection software samples data as it passes by the detector. Each “sampling” is stored as a data point. In ABI 373 and ABI PRISM 377 data, the scan number describes the location of the data point.

dye/sample Individual sample labeled with a single dye within a Sample file. Sample files normally contain up to four dye/samples, depending on how many labeled samples you included in each lane or injection of your data collection run.

dye color indicator Left color box in the Results Control window and the legend of the Results Display. In the Results Display, click this box to move the associated electropherogram to the front. In the Results Control and the Results Display windows, double-click this box to change the dye scale, or \mathbb{R} —double-click it to reset the dye scale to the default.

electropherogram Four-color picture of a sequence, showing peaks that represent the bases. The term is used interchangeably with chromatogram in this manual.

grid Spreadsheet-like display used for entering data in tabular format. The Analysis Control and Results Control windows display grids for entering sample information.

internal size standard Also called internal lane or injection size standard, DNA fragment of known size that you include with your run. On the ABI 373 or ABI PRISM 377, you include the size standard with the samples in each lane. On the ABI PRISM 310 you include the size standard with each injection.

Running an internal lane standard results in particularly accurate and precise molecular length determination because the internal lane standard and the unknown fragments undergo exactly the same electrophoretic forces. The software can then compensate for band-shift artifacts caused by variations in the gel and in the run. See “About Size Standards” on page 5-29 and “Using Size Standards” on page 5-36.

legend Informational text that appears beneath electropherogram panels in the results displays. You can show or hide legends (see “Using Legends to Change the Display” on page 7-30), and you can use the color boxes displayed in them to bring specified electropherograms to the front of the panel (“Moving the Electropherogram” on page 7-28) or customize the colors (“How to Define Custom Colors” on page 7-40).

matrix file/multicomponent matrix File used to adjust for the spectral overlap between the fluorescent dyes used on the ABI PRISM[®] instruments. A mathematical matrix of the spectral overlaps is created and the inverse matrix is used to correct the data during analysis. Matrix files are stored in the ABI folder inside the Macintosh[®] computer System folder, or in the matrix folder in the GeneScan Analysis Software folder. The values of the matrix are stored in the gel file (ABI 373 and ABI PRISM 377) and in the Sample files.

For more information, see Chapter 6, “Making a Matrix File.”

module A file that provides instructions about conditions of operation to the ABI PRISM genetic analysis instrument. Refer to your instrument manual.

multicomponenting Adjusting for spectral overlap of the fluorescent dyes.

overlaid Displayed together so they overlap. In the GeneScan Analysis Software Results Display window, all electropherograms in a single panel are overlaid. You can bring a specific one to the front by clicking the color box that represents it in the legend (See “Using Legends to Change the Display” on page 7-30).

plot color indicator Right color box in the Results Control window and the legend of the Results Display. In the Results Display window, click this box to move the associated electropherogram to the front. In the Results Control and the Results Display windows, double-click this box to change the plot color, or \mathbb{R} —double-click it to reset the plot color to the default.

preferences Defaults you can set so that certain parameters are automatically applied when you are working with a project. The GeneScan Analysis Software remembers preferences and applies them globally to all new projects.

For a brief description of each preference you can set, see “Changing How the Results are Displayed and Printed” on page 7-13.

project File containing links to a set of Sample files that you want to analyze and display together. A project can contain Sample files from multiple runs. Adding a Sample file to a project creates a reference to the file. It does not copy the file into the project.

For more information, see “Using a Project to Manage Sample Files” on page 4-8.

project options Formatting information you can set for the current project. Project options are remembered by the project when you open it again.

For a brief description of each project option, see “Changing How the Results are Displayed and Printed” on page 7-13.

sample files Computer files that contain raw and analyzed data. Sample files are created directly by the ABI PRISM 310 and by the GeneScan Analysis Software for the ABI 373 and ABI PRISM 377. Sample files contain data such as peak locations, size calling values, and a record of analysis settings.

scan number (ABI 373 and ABI PRISM 377 instruments) Data Collection software samples data 194 times (194 or 388 times for XL instruments) as it scans across the gel. Each “sampling” is stored as a data point. The scan number describes the location of the data point.

separation distance Length from the wells of the gel to the read region of the gel. Also called the “WTR (well-to-read).”

shark’s-tooth comb Piece of mylar inserted into a gel, typically used for a sequencing run. The flat edge of the comb is used to form a well during gel polymerization. The toothed edge is inserted to form wells into which samples are loaded. For GeneScan Analysis Software applications, a square-tooth comb is recommended to lessen the chance of sample spillage from lane to lane.

size calling curve Curve created by the GeneScan Analysis Software for size calling. The software calculates this curve based on the size calling method you specify for data analysis. This curve is red in the Standard Sizing Curve window. When it matches the align-by-size curve, the two overlap so you see only this curve. See also “align-by-size curve” and “size standard spline interpolation curve.”

size standard Specific DNA fragments of known sizes. After you define the peaks of a size standard, the GeneScan Analysis Software matches this definition to the internal lane or injection standard that you include with your run. The software assigns the defined size values to the appropriate peaks of the internal lane or injection standard, and uses this information with the selected size calling method to size all unknown fragments.

For more information, see “About Size Standards” on page 5-29 and “Using Size Standards” on page 5-36.

size standard spline interpolation curve Curve created by the GeneScan Analysis Software for aligning data by size. The software creates this curve if you use the Local Southern or Cubic Spline Interpolation size calling method and the size standard data does not match the best-fit curve, which is normally used for aligning the data by size. This curve is blue in the Sizing Curve window. See also “align-by-size curve” and “size calling curve.”

square-tooth comb Piece of mylar inserted into a gel to form wells for sample loading. You remove a square-tooth comb prior to sample loading. For GeneScan Analysis Software applications, a square-tooth comb is recommended to minimize the chance of lane-to-lane leakage, which can occur more often with a shark’s-tooth comb.

tiled Displayed so they do not overlap. The GeneScan Analysis Software displays tiled electropherogram panels in the Results Display. If you display more than one electropherogram in each panel, all electropherograms in the panel are overlaid.

tracker line Marks the channel exhibiting the strongest fluorescent signal in a lane. For each lane, the raw data extracted for the sample file is an average of the data in several channels. You set the number of channels averaged in the Gel Processing Parameters.

WTR (well-to-read) Length from the wells of the gel to the read region of the gel. See also “separation distance.”

Index

A

ABI 373

- automatic analysis diagram 4-3
- Collection file, opening A-7 to A-8
- gel processing parameters, setting A-5 to A-6
- generating sample files
 - procedure 6-27 to 6-28
 - when to generate new files A-12
- how GeneScan works with 1-5
- loading and running dye standards 6-20 to 6-23
- matrix standards, preparing 6-12
- processing using ABI 373 A-2 to A-4
- sample sheet, completing A-9 to A-10
- tracking lanes A-11

ABI 373 (XL), loading and running dye standards 6-24 to 6-26

ABI folder, files installed in 6-6

ABI PRISM 310

- automatic analysis diagram 4-2
- automatically printing run results 9-3
- loading and running dye standards 6-13 to 6-16
- matrix standards, preparing 6-9 to 6-11

ABI PRISM 377

- automatic analysis diagram 4-3
- automatically printing run results 9-3
- generating sample files 6-27 to 6-28
- loading and running dye standards 6-17 to 6-19
- matrix standards, preparing 6-12

Analysis Control window

- using to analyze project files 5-2 to 5-5
- using to analyze sample file 5-6 to 5-16

Analysis Log, using 7-51 to 7-52

analysis parameters

about 5-17 to 5-23

defined Glossary-1

using 5-24 to 5-28

analysis results

changing how displayed and printed 7-13 to 7-15

Results Display window

about 7-4 to 7-6

using 7-7 to 7-12

steps to evaluating

evaluating results 7-2

updating the results 7-16

using the Analysis Log 7-51 to 7-52

ways to display 7-3

See Also Sample Results View

analyzing

project files 5-2 to 5-5

sample files 5-6 to 5-16

archiving 8-6

automatic analysis, setting up 4-4 to 4-7

B

band-shift artifacts, compensating for Glossary-2

baselining defined Glossary-1

BioLIMS

about server names E-17 to E-19

accessing the database E-13 to E-16

configuring the server E-4 to E-9

GeneScan and BioLIMS, about E-2 to E-3

switching between Sample file and BioLIMS mode E-10 to E-11

testing network connection

using tool E-12

troubleshooting

client to Oracle connection F-14 to F-24

client to Sybase connection F-3 to F-13

GeneScan crashes with BioLIMS D-9
if BioLIMS Preference page does not
appear F-2
using the BioLIMS Chooser E-20 to E-30

C

channels defined Glossary-1
colors
 assigning to electropherograms 7-39
 defining in electropherograms 7-40 to 7-43
computer, using the Macintosh 1-7 to 1-10
configuring the BioLIMS server E-4 to E-9
copying data to other applications 8-7 to 8-8
crashes, GeneScan crashes with BioLIMS D-9
Cubic Spline method C-4
customer support
 e-mail address 1-23
 help 1-20 to 1-25
 internet address 1-20
 regional sales offices 1-23 to 1-25
 telephone/fax (U.S.) 1-20

D

Data Collection software
 automatically printing results from 9-3
 Gel File window, how differs from GeneScan
 Gel File window 2-12
data point defined Glossary-1
Deionizing formamide, procedure 6-9
dye color, changing in electropherograms 7-28
Dye Matrix Standards kits 6-8
dye scale, changing 7-44 to 7-45

E

electropherograms
 about 7-20 to 7-21
 changing dye scale 7-44 to 7-45
 defining custom colors 7-40 to 7-43
 displaying electropherogram data 7-25 to
 7-26
 displaying electropherograms and tabular
 data 7-22 to 7-24
 using to verify peak detection 7-53
 using to verify results 7-46 to 7-47
 verifying size calculations 7-48 to 7-50

working with electropherogram data 7-27 to
7-39
 assigning standard colors 7-39
 changing horizontal scale 7-37
 changing the dye color 7-28
 changing vertical scale 7-38 to 7-39
 for all electropherograms 7-38
 for single electropherograms 7-39
 displaying peak positions 7-29
 displaying x-and y-axis positions 7-27
 highlighting peaks 7-29
 making active window 7-28
 scrolling the display 7-30
 showing data by fragment size 7-35 to
 7-36
 about 7-35
 procedure 7-36
 showing off-scale data 7-32
 using legends to change the
 display 7-30
 zooming in and out 7-31
e-mail, address for technical support 1-23
EPT Data View 3-17 to 3-18
 displaying the view 3-17
 example 3-18
 what it displays 3-17
error messages D-9 to D-12
extracting data, and tracking 2-59 to 2-66
 extracting without changing tracking 2-64 to
 2-66
 saving after tracking 2-68
 specifying Run folder 2-59
 tracking and extracting 2-61 to 2-63
 tracking without extracting 2-60 to 2-61
 ways to track and extract 2-59

F

FAX on Demand 1-22
files, software files, table of 1-16 to 1-18
folders
 ABI folder, files installed in 6-6
 defining locations 5-41 to 5-43
formamide
 how to deionize, procedure 6-9
 size standard mix, preparing 6-10

G

- Gel File window
 - diagram 2-13
 - optimizing tracker lines 2-54
 - window callouts
 - call outs described 2-15 to 2-17
 - window buttons 2-14 to 2-15
 - gel files
 - about the gel files 2-2
 - archiving 8-6
 - displaying the gel file 2-11
 - Gel File window
 - diagram 2-13
 - differences between windows 2-12
 - window callouts
 - call outs described 2-15 to 2-17
 - window buttons 2-14 to 2-15
 - gel image, adjusting 2-18 to 2-25
 - adjusting the contrast 2-20 to 2-22
 - changing dye indicators 2-19 to 2-20
 - displaying/hiding a dye color 2-18
 - displaying/hiding dye color 2-18
 - installing new matrix file 2-24 to 2-25
 - regenerating the gel image 2-23 to 2-24
 - removing contrast changes 2-22
 - Gel Processing preferences, setting 2-3 to 2-10
 - Auto-Launch processing 2-4
 - comb type 2-10
 - displaying the Gel preferences 2-3
 - Image Generation Defaults 2-4 to 2-6
 - lane extraction 2-7 to 2-10
 - interpolating data 2-57 to 2-58
 - lane markers, working with 2-43 to 2-52
 - off-scale data, displaying 2-38 to 2-41
 - sample sheet, about 2-29 to 2-37
 - saving
 - after editing tracking 2-68
 - procedure 8-4
 - tracker lines, working with 2-53 to 2-56
 - tracking and extracting data 2-59 to 2-66
 - extracting without changing tracking 2-64 to 2-66
 - saving after tracking 2-68
 - specifying Run folder 2-59
 - tracking and extracting 2-61 to 2-63
 - tracking without extracting 2-60 to 2-61
 - ways to track and extract 2-59
 - troubleshooting gel data D-5 to D-6
 - verifying gel information 2-26 to 2-28
 - why save 8-2
- gel image
 - adjusting 2-18 to 2-25
 - adjusting the contrast 2-20 to 2-22
 - changing dye indicators 2-19 to 2-20
 - displaying/hiding a dye color 2-18
 - displaying/hiding gel color 2-18
 - regenerating the gel image 2-23 to 2-24
 - removing contrast changes 2-22
 - printing 9-4
 - GeneScan
 - about this manual 1-5 to 1-6
 - archiving 8-6
 - copying, data to other applications 8-7 to 8-8
 - crashes with BioLIMS D-9
 - license and warranty G-1 to G-3
 - preparing matrix standard samples 6-10
 - registering 1-11
 - saving
 - gel files 8-4
 - projects 8-3
 - Results Displays 8-5
 - sample files 8-3
 - why save 8-2
 - size standard
 - GeneScan 1000 B-8 to B-9
 - GeneScan 2500 B-10 to B-11
 - GeneScan 350 B-2 to B-3
 - GeneScan 400HD B-4 to B-5
 - GeneScan 500 B-6 to B-7
 - software files, table of 1-16 to 1-18
 - what new in this release 1-2 to 1-3
 - GeneScan 100 standard B-8 to B-9
 - GeneScan 2500 standard B-10 to B-11
 - GeneScan 350 size standard B-2 to B-3
 - GeneScan 500 size standard B-6 to B-7
 - GeneScan-400 HD size standard B-4 to B-5
 - Genotyper
 - copying GeneScan data to 8-7
 - troubleshooting software results D-8
 - Global Southern method C-7 to C-8
 - glossary Glossary-1 to Glossary-4
 - grid defined Glossary-2

H

help

- e-mail address 1-23
 - internet address 1-20
 - regional sales offices 1-23 to 1-25
 - telephone hours 1-20
 - telephone/fax (U.S.) 1-20
- horizontal scale, changing in
- electropherograms 7-37

I

- IMPORTANT, definition 1-6
- installing 1-12 to 1-15
- folder contents 1-15
 - procedure 1-14
 - program files installed, diagram 1-15
 - RAM required 1-14
 - system specifications 1-12
- instrument, manual text related to 1-5
- internal size standard
- compensating for band-shift artifacts Glossary-2
 - defined Glossary-2
- internet address
- FAX-on-Demand 1-22
 - Perkin-Elmer 1-20
- interpolating data 2-57 to 2-58

L

- lane markers, working with 2-43 to 2-52
- Least Square size calling method C-2 to C-3
- legends
- defined Glossary-2
 - using to change display in electropherograms 7-30
- license and warranty G-1 to G-3
- Local Southern method C-5 to C-6

M

- Macintosh, using the computer 1-7 to 1-10
- manual, text for specific instruments 1-5
- matrix files
- about 6-2 to 6-6
 - assigning to sample files 6-35 to 6-36
 - bad matrix files, causes 6-38 to 6-39
 - choosing a scan range 6-29 to 6-31

- defined Glossary-2
 - Dye Matrix Standard kits 6-8
 - evaluating the matrix file 6-37
 - generating files (ABI 373 and ABI PRISM 377) 6-27 to 6-28
 - installing new file
 - gel file 2-24 to 2-25
 - sample file 3-20
 - loading and running dye standards
 - ABI 373 6-20 to 6-23
 - ABI 373 (XL) 6-24 to 6-26
 - ABI PRISM 310 6-13 to 6-16
 - ABI PRISM 377 6-17 to 6-19
 - new matrix file, generating 6-32 to 6-33
 - preparing matrix standards
 - ABI PRISM 310 6-9 to 6-11
 - ABI Prism 377 and ABI 373 6-12
 - process of creating new file 6-7 to 6-8
 - saving and naming 6-34
- matrix standards, preparing standard samples
- GeneScan 6-10
- module defined Glossary-2
- multicomponent matrix defined Glossary-2
- multicomponenting defined Glossary-2

N

- network, using SybPing program E-12
- note, definition 1-6

O

- off-scale data
- showing in electropherograms 7-32
 - showing in gel files 2-38 to 2-41
- optimizing tracker lines 2-54
- Oracle server
- configuring the server E-4 to E-9
 - how server names are recognized E-17 to E-19
 - troubleshooting connection F-14 to F-24
- overlaid defined Glossary-2

P

- PCR Single-stranded Conformation Polymorphism (SSCP)
- Minimum Peak Half width setting 5-21
- peak detection, verifying 7-53

- peak positions, displaying in
 - electropherograms 7-29
 - peaks, highlighting in electropherograms 7-29
 - preferences, defined Glossary-3
 - printing
 - about printing 9-2
 - automatically printing run results 9-3
 - gel image 9-4
 - sample files 9-5 to 9-6
 - supporting information 9-6
 - program files, installed (diagram) 1-15
 - project options, defined Glossary-3
 - projects
 - analyzing project files 5-2 to 5-5
 - creating
 - finding missing sample files 4-14 to 4-15
 - process diagrams 4-2 to 4-3
 - using ABI 373 and ABI PRISM 377 4-3
 - using the ABI PRISM 310 4-2
 - using to manage sample files 4-8 to 4-9
 - working with 4-10 to 4-13
 - creating a new project 4-10 to 4-12
 - opening an existing project 4-10
 - removing samples from a project 4-13
 - defined Glossary-3
 - defining folder locations 5-41 to 5-43
 - saving, procedure 8-3
 - troubleshooting D-2 to D-4
 - why save 8-2
- R**
- RAM required 1-14
 - Raw Data View 3-15 to 3-16
 - displaying the view 3-15
 - example 3-16
 - what it displays 3-15
 - what to evaluate 3-16
 - registering the software 1-11
 - restarting Macintosh, key commands for E-12
 - Results Control window
 - saving and renaming 7-17 to 7-19
 - important considerations 7-17
 - renaming current display 7-19
 - saving the format 7-17
 - using previously saved formats 7-18
 - working with previously saved displays 7-18
 - Results Display window
 - about 7-4 to 7-6
 - using 7-7 to 7-12
 - Results Displays
 - saving, procedure 8-5
 - why save 8-2
- S**
- Sample File window
 - about 3-5
 - EPT Data View 3-17 to 3-18
 - displaying the view 3-17
 - example 3-18
 - what it displays 3-17
 - Raw Data View 3-15 to 3-16
 - displaying the view 3-15
 - example 3-16
 - what it displays 3-15
 - what to evaluate 3-16
 - Sample Info View 3-8 to 3-12
 - description of view 3-9 to 3-12
 - displaying the view 3-8
 - example 3-9
 - what it displays 3-8
 - Sample Results View 3-6 to 3-7
 - description of view 3-7
 - differences from Results Display 3-7
 - displaying the view 3-6
 - example 3-6
 - what it displays 3-6
 - Size Curve View 3-13 to 3-14
 - displaying the view 3-13
 - size curve described 3-14
 - what is the size curve 3-13
 - what the size curve displays 3-14
 - sample files
 - about 3-2 to 3-3
 - how GeneScan analyzes sample files 3-3
 - how they are generated 3-2
 - ways to generate sample files 3-3
 - what sample files contain 3-2
 - analyzing 3-19 to 3-20
 - installing new matrix file 3-20
 - procedure 3-19 to 3-20

- using Analysis Control window 5-6 to 5-16
- archiving 8-6
- defined Glossary-3
- finding missing sample files 4-14 to 4-15
- how GeneScan names files 2-67
- matrix file, assigning to sample files 6-35 to 6-36
- opening, procedure 3-4
- printing 9-5 to 9-6
- removing from a project 4-13
- Sample File window
 - about 3-5
 - EPT Data view 3-17 to 3-18
 - Raw Data View 3-15 to 3-16
 - Sample Info View 3-8 to 3-12
 - Sample Results View 3-6 to 3-7
 - Size Curve View 3-13 to 3-14
- saving, procedure 8-3
- switching between Sample file and BioLIMS mode E-10 to E-11
- unlocking sample files 4-12
- using project to manage files 4-8 to 4-9
- why save 8-2
- Sample Info View 3-8 to 3-12
 - description of view 3-9 to 3-12
 - displaying the view 3-8
 - example 3-9
 - what it displays 3-8
- Sample Results View 3-6 to 3-7
 - description of view 3-7
 - differences from Results Display 3-7
 - displaying the view 3-6
 - example 3-6
 - what it displays 3-6
- sample sheet
 - about 2-29 to 2-37
 - completing for the ABI 373 A-9 to A-10
- saving
 - after tracking 2-68
 - gel files 8-4
 - matrix file 6-34
 - projects 8-3
 - Results Control format 7-17 to 7-19
 - important considerations 7-17
 - renaming current display 7-19
 - renaming current Results Display 7-19
 - saving 7-17
 - saving the format 7-17
 - using previously saved formats 7-18
 - working with previously saved displays 7-18
 - Results Displays 8-5
 - sample files 8-3
 - why save 8-2
- scan numbers, defined Glossary-3
- separation distance, defined Glossary-3
- server names, about E-17 to E-19
- sharks-tooth comb, defined Glossary-3
- size calling methods
 - Cubic Spine method C-4
 - Global Southern method C-7 to C-8
 - Least Square method C-2 to C-3
 - Local Southern method C-5 to C-6
- Size Curve View 3-13 to 3-14
 - displaying the view 3-13
 - size curve described 3-14
 - what is the size curve 3-13
 - what the size curve displays 3-14
- size standard
 - about 5-29 to 5-30
 - defining, how to 5-31 to 5-35
 - GeneScan 1000 B-8 to B-9
 - GeneScan 2500 B-10 to B-11
 - GeneScan 350 B-2 to B-3
 - GeneScan 400HD B-4 to B-5
 - GeneScan 500 B-6 to B-7
 - using 5-36 to 5-40
- Slice view, how to determine the scale of the peaks in view 2-6
- software
 - files, table of 1-16 to 1-18
 - license G-1 to G-3
 - registering 1-11
- SSCP, Minimum Peak Half width setting 5-21
- Sybase SQL server
 - configuring the BioLIMS server E-4 to E-9
 - how server names are recognized E-17 to E-19
 - troubleshooting connection F-3 to F-13
- SybPing, using E-12
- system specifications 1-12

T

- tabular data displays
 - about 7-20 to 7-21
 - displaying electropherogram and tabular data 7-22 to 7-24
 - using to verify peak detection 7-53
 - using to verify results 7-46 to 7-47
 - technical support 1-20 to 1-25
 - e-mail address 1-23
 - internet address 1-20
 - regional sales offices 1-23 to 1-25
 - telephone/fax (U.S.) 1-20
 - text in manual, noted for specific instruments 1-5
 - tracking, and extracting data
 - extracting without changing tracking 2-64 to 2-66
 - saving after tracking 2-68
 - specifying Run folder 2-59
 - tracker lines
 - defined Glossary-4
 - working with 2-53 to 2-56
 - tracking and extracting 2-61 to 2-63
 - tracking without extracting 2-60 to 2-61
 - ways to track and extract 2-59
- troubleshooting
- bad matrix files, causes 6-38 to 6-39
 - BioLIMS
 - client to Oracle connection F-14 to F-24
 - client to Sybase connection F-3 to F-13
 - GeneScan crashes with BioLIMS D-9
 - if BioLIMS Preference page does not appear F-2
 - error messages D-9 to D-12
 - gel data D-5 to D-6
 - Genotyping software results D-8
 - projects and results D-2 to D-4

U

- unlocking sample files 4-12

V

- vertical scale, changing 7-38 to 7-39
 - for all electropherograms 7-38
 - for single electropherograms 7-39

- viewing scale, changing for electropherograms 7-31

W

- warranty G-1 to G-3
- well-to-read (WTR), defined Glossary-4
- www address
 - FAX-on-Demand 1-22
 - Perkin-Elmer 1-20

X

- x-axis positions, displaying in electropherograms 7-27

Y

- y-axis positions, displaying in electropherograms 7-27

Z

- zooming, electropherograms 7-31

