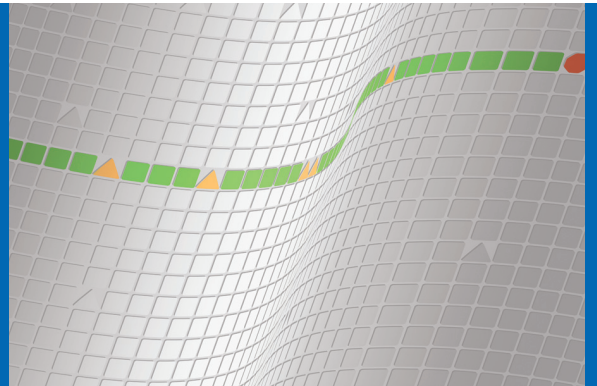


# GeneMapper® *ID-X* Software

## Version 1.2

**Note:** To improve the clarity of graphics in this PDF file, use the zoom tool to increase magnification to 150% or greater.

GeneMapper®  
***ID-X***



© Copyright 2009, Life Technologies Corporation. All rights reserved.

**For Research, Forensic, or Paternity Use Only. Not for use in diagnostic procedures.**

Information in this document is subject to change without notice. Life Technologies Corporation assumes no responsibility for any errors that may appear in this document.

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

**Notice to Purchaser: License Disclaimer.**

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

GeneMapper® *ID-X* Software has undergone a verification process defined by Applied Biosystems. However, human identification laboratories analyzing forensic, paternity, databasing, and single-source samples that choose to use GeneMapper® *ID-X* Software for data analysis should perform their own appropriate validation studies.

**TRADEMARKS:**

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

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

Part Number 4426481 Rev. A  
06/2009

# Contents

Preface	vii
Revision History	vii
How to Use This Guide	vii
Purpose of This Guide	vii
Audience	vii
Assumptions	vii
Text Conventions	vii
User Attention Words	viii
Send Us Your Comments	viii
How to Obtain Support	ix
Chapter 1	
Peak Detection, Sizing, Allelic Ladder, and Sample Analysis Algorithms	1
Allelic Ladder Requirements	2
Allelic Ladder Run Requirements	2
Allelic Ladder Analysis Requirements	3
Overview of the Peak Detection, Sizing, and Genotyping Process	4
Peak Detection	4
Examining Peak Definitions	5
Baseline Window Size	6
Guideline for Using This Parameter	6
Effects of Varying Baseline Window Size	6
Smoothing	8
Polynomial Degree and Peak Window Size	9
Effects of Varying the Polynomial Degree	9
Effects of Increasing the Window Size Value	10
Optimizing Peak Detection Sensitivity – Example 1: Reducing Peak Window Size	11
Effects of Reducing Window Size	12

Optimizing Peak Detection Sensitivity – Example 2: Extreme Settings .....	13
Slope Thresholds for Peak Start/End Parameters .....	13
Guidelines for Using These Parameters .....	14
Using These Parameters .....	14
Slope Threshold Example – Asymmetrical Peak .....	15
Adjusting Slope Threshold for Peak End .....	15
Sizing .....	16
Size Matching .....	16
Size-Calling Curve Generation and Size Calling .....	18
Size Calling – Local Southern Method .....	19
Local Southern Method Equation .....	19
How The Local Southern Method Works .....	20
Size Calling – Global Southern Method .....	21
Global Southern Method Equations .....	21
How the Global Southern Method Works .....	21
Size Calling – Least Squares Method .....	22
Using 3rd Order for MiniFiler™ Kit Data .....	22
Examples .....	22
Size Calling – Cubic Spline Interpolation Method .....	24
Possible Local Sizing Inaccuracy .....	25
Sizing Quality Assessment .....	25
Allelic Ladder Genotyping .....	26
Allele Matching .....	26
Allelic Ladder Quality Assessment .....	27
Bin Offsetting .....	27
How the Software Calculates and Applies Bin Offsets .....	29
Example: One Allelic Ladder Is Present In a Run Folder .....	30
Example: More Than One Allelic Ladder Sample Is Present In a Run Folder .....	32
Example: How Offsets Are Applied To Virtual Alleles .....	32
Displaying Offset Marker Margins .....	33

---

	Sample Genotyping	34
	Allele Calling and Filtering of Samples	34
	Off Ladder (OL) Allele Calls	35
	Allele Naming Convention	35
	Genotype and Sample Quality Assessment	36
<b>Chapter 2</b>	<b>GeneMapper® <i>ID-X</i> Software Quality Value System</b>	<b>37</b>
	Overview of the Quality Value System	38
	Optimizing and Validating an Expert System	38
	Displaying Numeric SQ and GQ Values	39
	Overview of the Quality Value System Checks and Assessments	40
	Analysis Requirement Checks	41
	Sizing Quality Assessment	43
	Allelic Ladder Quality Assessment	45
	Quality Rules	45
	Effect on GQ	46
	Marker-Level Quality Assessment	47
	Genotype Quality Assessment	51
	Sample-Level Quality Assessments	53
<b>Appendix A</b>	<b>Troubleshooting</b>	<b>57</b>
	Software Issues	57
	<b>Bibliography</b>	<b>59</b>
	<b>Index</b>	<b>61</b>



# Preface

## Revision History

Part Number	Revision	Date	Description of Changes
4426481	A		New Document for Version 1.2 Software: Improvements to allelic ladder algorithm

## How to Use This Guide

**Purpose of This Guide** The *GeneMapper*<sup>®</sup> *ID-X* Software Version 1.2 Reference Guide:

- Provides reference information on sizing and genotyping.
- Describes the *GeneMapper*<sup>®</sup> *ID-X* Software quality value system and peak quality values (PQVs).

**Audience** This guide is intended for *GeneMapper ID-X* Software users who review and evaluate sizing and genotyping data.

**Assumptions** This guide assumes that you are familiar with analyzing data using the *GeneMapper*<sup>®</sup> *ID-X* Software.

**Text Conventions** This guide uses the following conventions:

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

- A right arrow bracket ( ▶ ) separates successive commands you select from a drop-down or shortcut menu. For example:  
Select **File ▶ Open ▶ Spot Set**.  
Right-click the sample row, then select **View Filter ▶ View All**.

## User Attention Words

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

---

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

---

---

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

---

Examples of the user attention words appear below:

---

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

---

---

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

---

---

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

---

---

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

---

## Send Us Your Comments

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

[techpubs@appliedbiosystems.com](mailto:techpubs@appliedbiosystems.com)



## How to Obtain Support

For HID support, you can send an e-mail to [HIDTechSupport@appliedbiosystems.com](mailto:HIDTechSupport@appliedbiosystems.com) or call **888-821-4443** option **1**.

For HID support outside North America, contact your local support office.

For the latest services and support information for all locations, go to <http://www.appliedbiosystems.com>, then click **Support**. At the Support page, you can:

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



## 1

# Peak Detection, Sizing, Allelic Ladder, and Sample Analysis Algorithms

1

This chapter covers:

- Allelic Ladder Requirements. . . . .2
- Overview of the Peak Detection, Sizing, and Genotyping Process. . . . .4
- Peak Detection. . . . .4
- Sizing. . . . .16
- Allelic Ladder Genotyping . . . . .26
- Sample Genotyping . . . . .34

## Allelic Ladder Requirements

### Allelic Ladder Run Requirements

To accurately genotype samples, you must use an Allelic Ladder sample that has been run with the samples. For samples that are run on an:

- ABI PRISM<sup>®</sup> 310 Genetic Analyzer – Run at least one allelic ladder for every 10 sample injections.
- ABI PRISM<sup>®</sup> 3100 or Applied Biosystems 3130 series Genetic Analyzers – Run at least one allelic ladder for **every set of 16 samples**.
  - 3130*x*l or 3100 Genetic Analyzers – One ladder per injection (one injection = 16 samples)
  - 3130 or 3100-Avant Genetic Analyzers – One ladder for every four injections (one injection = 4 samples)

---

**IMPORTANT!** Variation in laboratory temperature can cause changes in fragment migration speed that can, in turn, cause sizing variation. Applied Biosystems recommends the frequency of allelic ladder injections described above to account for normal variation in fragment migration speed. However, during internal validation studies, verify the required allelic ladder injection frequency to ensure accurate genotyping of all samples in your laboratory environment.

---

It is critical to genotype using an allelic ladder that is run under the same conditions as the samples because:

- Size values that are obtained for the same sample can differ between instrument platforms because of different polymer matrices and electrophoretic conditions.
- Slight procedural and reagent variations between single and multiple capillaries result in greater size variation than that found between samples injected in the same capillary in a single run.

## Allelic Ladder Analysis Requirements

- To accurately genotype samples, the GeneMapper® *ID-X* Software requires at least one allelic ladder sample per run folder.

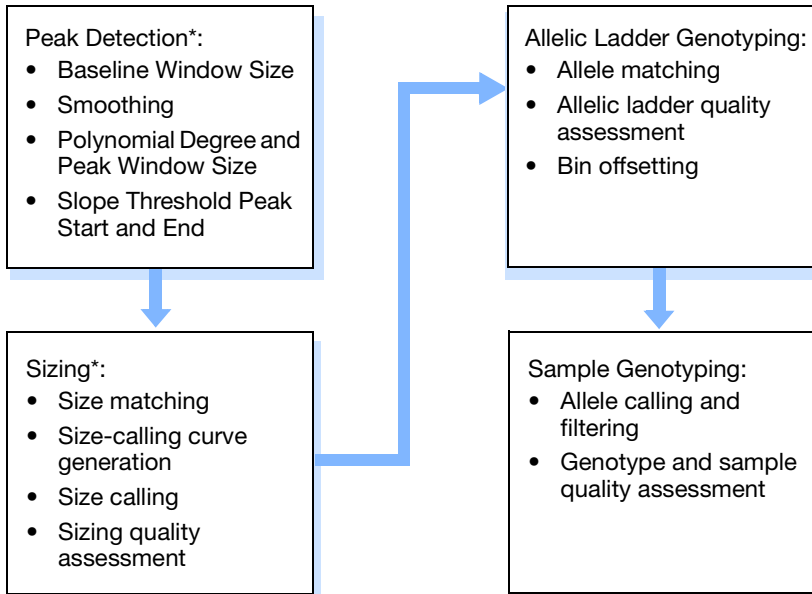
---

**IMPORTANT!** If a run folder contains multiple allelic ladder samples, the GeneMapper *ID-X* Software calculates bin offsets using an average of all ladders that use the same panel within a run folder (for more information, see [“Bin Offsetting” on page 27](#)).

---

- Allelic ladder samples must be identified as “Allelic Ladder” in the Sample Type column in a project.
- Analyze the allelic ladder sample(s) using the same analysis method and parameters used for samples.

## Overview of the Peak Detection, Sizing, and Genotyping Process



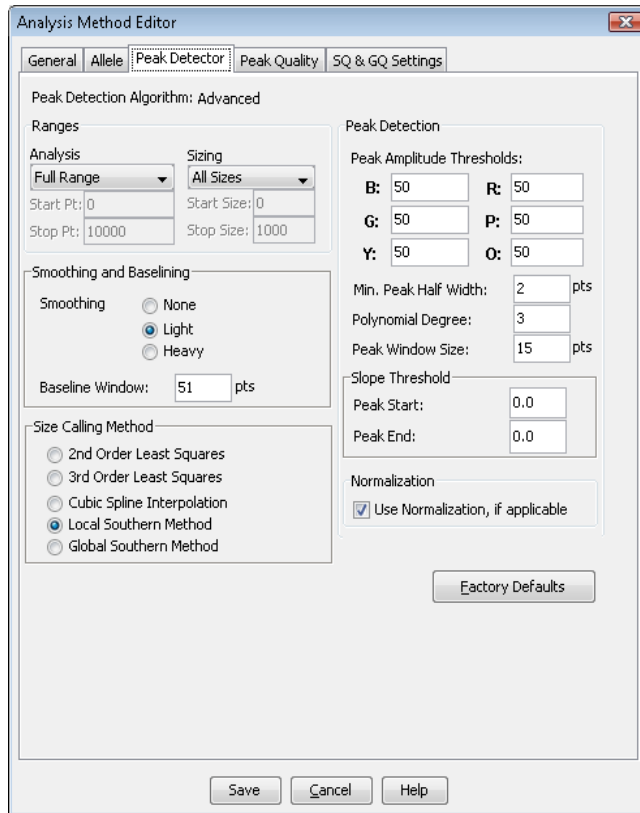
\*Sample peak detection and sizing is performed after allelic ladder genotyping.

## Peak Detection

### Peak Detection:

- Baseline Window Size
- Smoothing
- Polynomial Degree and Peak Window Size
- Slope Threshold Peak Start and Peak End

Parameters in the Analysis Method Peak Detector tab determine how the raw data are baselined and smoothed, and how peaks are detected.



This section describes how these parameters work. For more information on these parameters, see the *GeneMapper® ID-X Software Help*.

**IMPORTANT!** This section provides examples of the effect of changing certain peak detection parameters. Applied Biosystems provides default settings in the software, but your laboratory should verify these settings, or changes to these settings, as part of your laboratory’s internal validation studies.

## Examining Peak Definitions

To examine how GeneMapper® ID-X Software has defined a peak, select **Plot ▶ Peak Positions** in the Samples plot window. The peak positions (start, apex, end) are tick-marked in the electropherogram.

## Baseline Window Size

### Peak Detection:

- **Baseline Window Size**
- Smoothing
- Polynomial Degree and Peak Window Size
- Slope Threshold Peak Start and Peak End

The Baseline Window size parameter controls baselining for a group of peaks.

The software determines a reference baseline value for each data point. In general, the software sets the reference baseline to be the lowest value that it detects in a specified window size (in data points) centered on each data point.

A small baseline window relative to the width of a cluster, or grouping of peaks spatially close to each other, can result in shorter peak heights.

Larger baseline windows relative to the peaks being detected can create an elevated baseline, resulting in peaks that are elevated or not resolved to the baseline.

### Guideline for Using This Parameter

The default setting of 51 is optimal for analysis of AmpF $\Phi$ STR<sup>®</sup> kit data.

Choose a value that encompasses the width in data points of the peaks being detected while preserving a qualitatively smooth baseline.

The trade-off for a smoother baseline that touches all peaks is a reduction in peak height.

### Effects of Varying Baseline Window Size

Figure 1 on page 7 shows a sample with different reference baselines (zero in the analyzed electropherogram) that result from different baseline window size settings:

- The red trace shows a baseline derived from an extreme baseline window size value of 2801. At this setting, the reference baseline does not touch all peaks and elevates peak heights.
- The blue trace shows a baseline derived from the default value of 51 data points.



- The black trace shows a baseline that is derived from an extreme baseline window size value of 5 data points. At this setting, the reference baseline tracks the peaks, significantly reducing peak height.

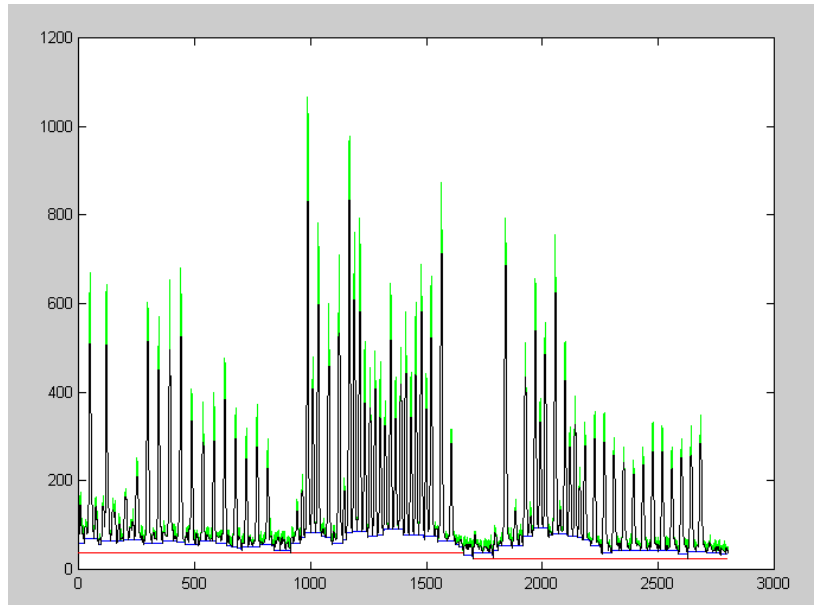


Figure 1 Baseline Window example

## Smoothing

### Peak Detection:

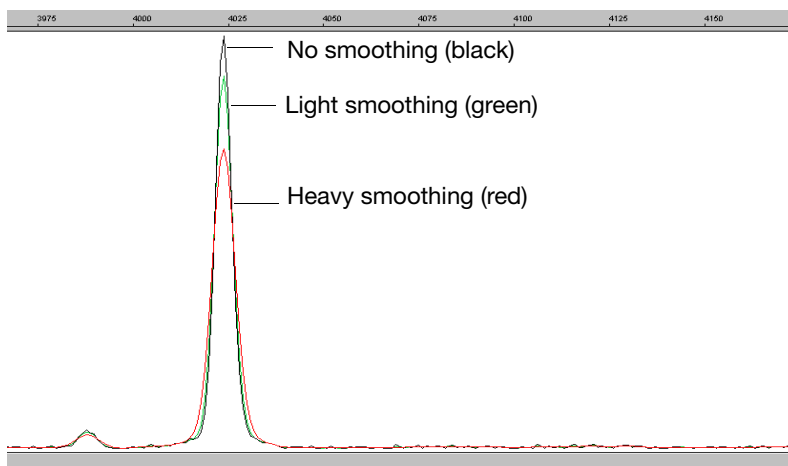
- Baseline Window Size
- **Smoothing**
- Polynomial Degree and Peak Window Size
- Slope Threshold Peak Start and Peak End

This parameter smooths the outline of peaks, and reduces the number of false peaks that are detected.

Smoothing is performed before peak detection and can be set to:

Option	Description
None	Applies no smoothing. Select for slower runs with very broad peaks, or to avoid the detection of sharp edges.
Light	Provides the best results for typical data. Light smoothing slightly reduces peak height.
Heavy	Select for data with very sharp, narrow peaks of interest. Heavy smoothing can significantly reduce peak height.

Figure 2 shows the effect of smoothing settings on a peak.



**Figure 2** Smoothing example

## Polynomial Degree and Peak Window Size

- Peak Detection:
- Baseline Window Size
  - Smoothing
  - **Polynomial Degree and Peak Window Size**
  - Slope Threshold Peak Start and Peak End

Polynomial Degree and Peak Window Size settings affect the peak detection sensitivity. You can adjust these parameters to detect a single base pair difference while minimizing the detection of shoulder effects and/or noise.

The peak window size functions with the polynomial degree to set the sensitivity of peak detection. The peak detector calculates the first derivative of a polynomial curve fitted to the data within a window that is centered on each data point in the analysis range.

Using curves with larger polynomial degree values allows the curve to more closely approximate the signal and, therefore, the peak detector captures more of the peak structure in the electropherogram.

The peak window size sets the width (in data points) of the window to which the polynomial curve is fitted to data:

- Higher peak window size values smooth out the polynomial curve, which limits the structure being detected.
- Smaller window size values capture more of the peak structure.

### How to Use the Peak Detection Parameters

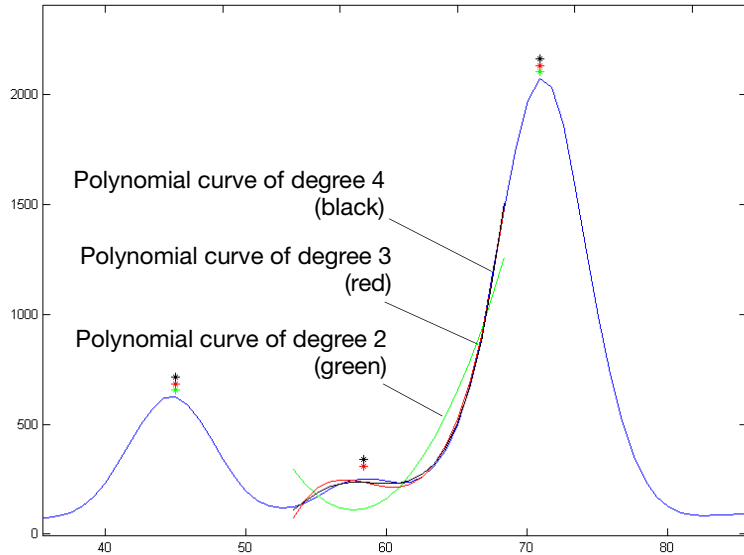
Use the table below to adjust the sensitivity of detection.

Function	Polynomial Degree Value	Window Size Value
Increase sensitivity	Higher	Lower
Decrease sensitivity	Lower	Higher

### Effects of Varying the Polynomial Degree

Figure 3 on page 10 shows peaks detected with a window size of 15 data points and a polynomial curve of degree 2 (green), 3 (red), and 4 (black). The diamonds represent a detected peak using the respective polynomial curves.

Note that the smaller trailing peak is not detected using a degree of 2 (green). As the peak detection window is applied to each data point across the displayed region, a polynomial curve of degree 2 could not be fitted to the underlying data to detect its structure.



**Figure 3** Electropherogram showing peaks detected with three different polynomial degrees

### Effects of Increasing the Window Size Value

In [Figure 4 on page 11](#) both polynomial curves have a degree of 3 and the window size value was increased from 15 (red) to 31 (black) data points.

As the cubic polynomial is stretched to fit the data in the larger window size, the polynomial curve becomes smoother. Note that the structure of the smaller trailing peak is no longer detected as a distinct peak from the adjacent larger peak to the right.

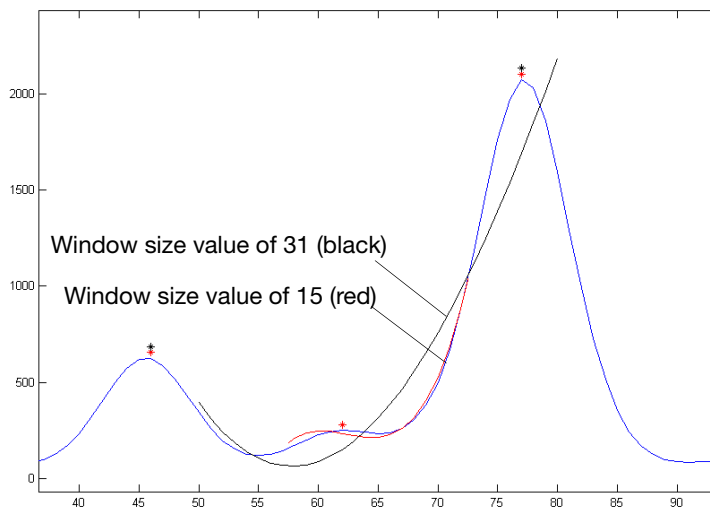


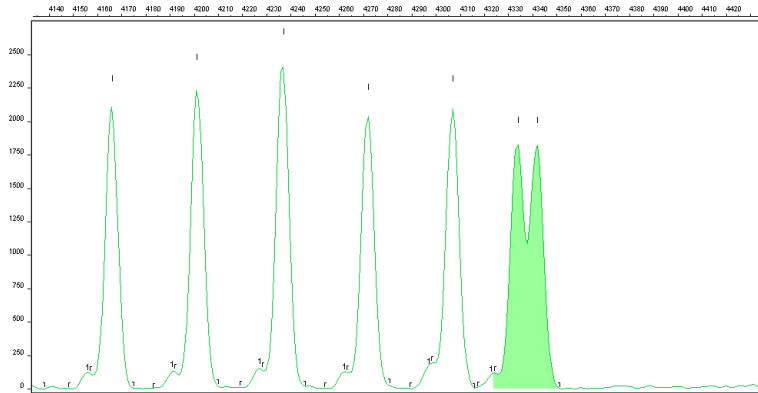
Figure 4 Electropherogram showing the same peaks as in Figure 3 after increasing the window size value but keeping the polynomial degree the same

## Optimizing Peak Detection Sensitivity – Example 1: Reducing Peak Window Size

### Peak Detection:

- Baseline Window Size
- Smoothing
- **Polynomial Degree and Peak Window Size**
- Slope Threshold
- Peak Start and Peak End

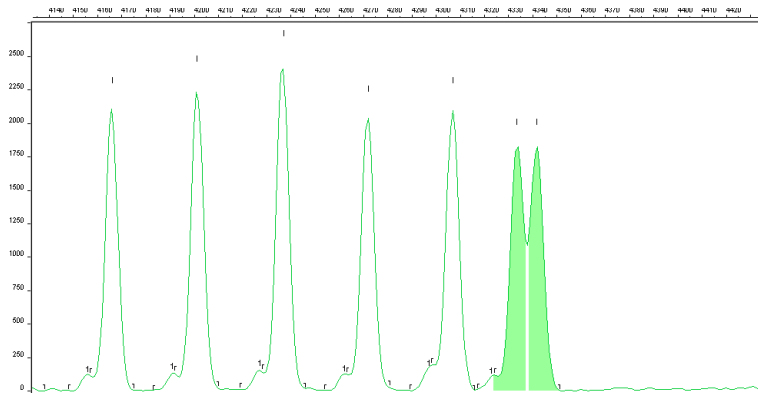
Figure 5 on page 12 shows the initial electropherogram with two resolved alleles of known fragment lengths (they differ by one nucleotide) that are detected as a single peak. The analysis was performed using a polynomial degree of 3 and a peak window size of 19 data points.



**Figure 5** Electropherogram showing two resolved alleles detected as a single peak

### Effects of Reducing Window Size

**Figure 6** shows that both alleles are detected after reanalyzing with the polynomial degree set to 3 and the window size value decreased to 15 (from 19) data points.



**Figure 6** Electropherogram showing the alleles detected as two peaks after decreasing the window size value

## Optimizing Peak Detection Sensitivity – Example 2: Extreme Settings

- Peak Detection:
- Baseline Window Size
  - Smoothing
  - **Polynomial Degree and Peak Window Size**
  - Slope Threshold Peak Start and Peak End

Figure 7 shows the result of an analysis using a peak window size value set to 10 and a polynomial degree set to 9. These extreme settings for peak detection caused several peaks to be split and detected as two separate peaks.

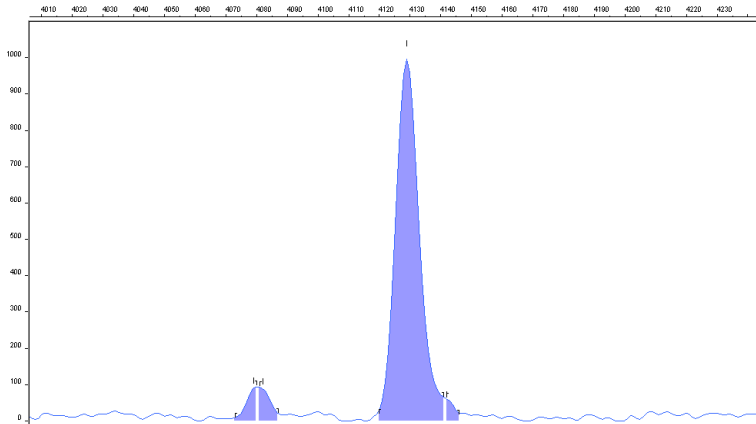


Figure 7 Electropherogram showing the result of an analysis using extreme setting for peak detection

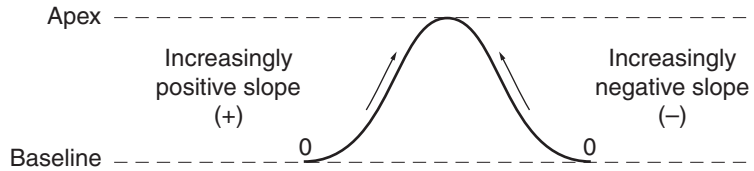
## Slope Thresholds for Peak Start/End Parameters

- Peak Detection:
- Baseline Window Size
  - Smoothing
  - Polynomial Degree and Peak Window Size
  - **Slope Threshold Peak Start and Peak End**

The Slope Threshold for Peak Start and Slope Threshold for Peak End parameters adjust the start and end points of a peak.

The values assigned to these parameters can be used to better position the start and end points of an asymmetrical peak, or a poorly resolved shouldering peak to more accurately reflect the peak position and area.

In general, from left to right, the slope of a peak increases from the baseline up to the apex. From the apex down to the baseline, the slope decreases negatively until it returns to zero at the baseline (Figure 8 on page 14).



**Figure 8** Peak slope

If either of the slope values you enter exceeds the slope of the peak being detected, the software overrides your value and reverts to zero.

### Guidelines for Using These Parameters

- For typical or symmetrical peaks, use a value of zero.
- For asymmetrical peaks, select values other than zero to better reflect the beginning and end points.
- A value of zero does not affect the sizing accuracy or precision of an asymmetrical peak.

### Using These Parameters

**Note:** The size of a detected peak is the calculated apex between the start and end points of a peak. Peak size does not change based on start and end settings.

To move the...	Then...	Example
Start point of a peak closer to its apex	Change the Slope Threshold for Peak Start value from zero to a positive number.	
End point of a peak closer to its apex	Change the Slope Threshold for Peak End value to a more negative number.	



## Slope Threshold Example – Asymmetrical Peak

- Peak Detection:
- Baseline Window
  - Smoothing
  - Polynomial Degree and Peak Window Size
  - **Slope Threshold Peak Start and Peak End**

### Initial Electropherogram

Figure 9 shows the initial electropherogram analyzed with value of 0 for Peak Start and Peak End. Note the asymmetrical peak with a noticeable tail on the right side.

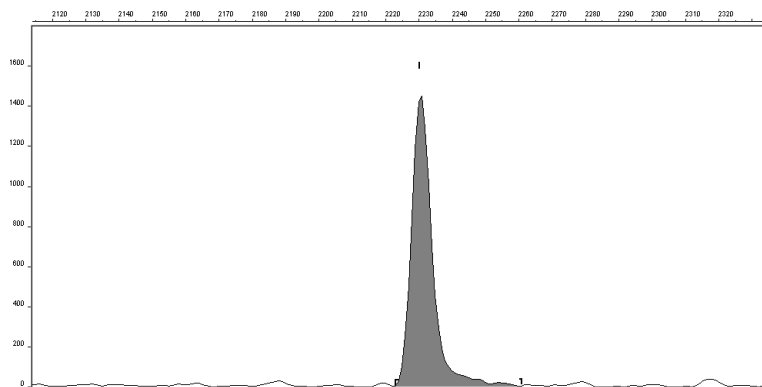


Figure 9 Electropherogram showing an asymmetrical peak

### Adjusting Slope Threshold for Peak End

After reanalysis with a value of  $-35.0$  for the Slope Threshold for Peak End, the end point that defines the peak moves closer to its apex, thereby removing the tail (Figure 10). Note that the only change to tabular data is the area (peak size and height are unchanged).

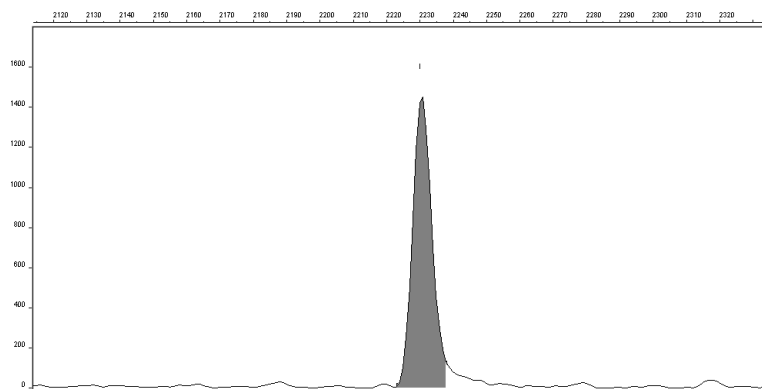


Figure 10 Electropherogram showing the effect of changing the slope threshold for peak end

## Sizing

**Sizing:**

- Size matching
- Size-calling curve generation
- Size calling
- Sizing quality assessment

During sizing, the software:

- Performs size matching of the internal size standard in all samples against the size standard definition selected in the software.
- Generates a size-calling curve.
- Sizes DNA sample peaks.
- Assesses the sizing quality.

## Size Matching

**Sizing:**

- **Size matching**
- Size-calling curve generation
- Size calling
- Sizing quality assessment

During size matching, the software matches the size standard fragments from the electropherogram to the list of fragment sizes in the size standard definition specified in the software.

Size matching uses ratio matching, based on relative height and distance of neighboring peaks. It then derives quality values statistically by examining the similarity between the theoretical (from the size standard definition) and actual (observed) fragment patterns.

The software ignores anomalous peaks that do not match the expected patterns. The software constructs a best-fit curve using the data points of each size standard fragment detected. A comparison between the sizes calculated from the best-fit curve and the matched peaks from the size standard definition file using the array of numbers is performed. Size-matching (and subsequent size calling) fails if significant differences in peak patterns are found, if no match can be made based on the expected patterns, or if all peaks are not found.

Because the software uses ratio-matching (looks for the expected number of alleles and expected peak patterns instead of specific data points), it is not necessary to define new size-standard definitions due to migration shifts.

[Figure 11 on page 17](#) illustrates the size matching and size-calling curve generation algorithm.

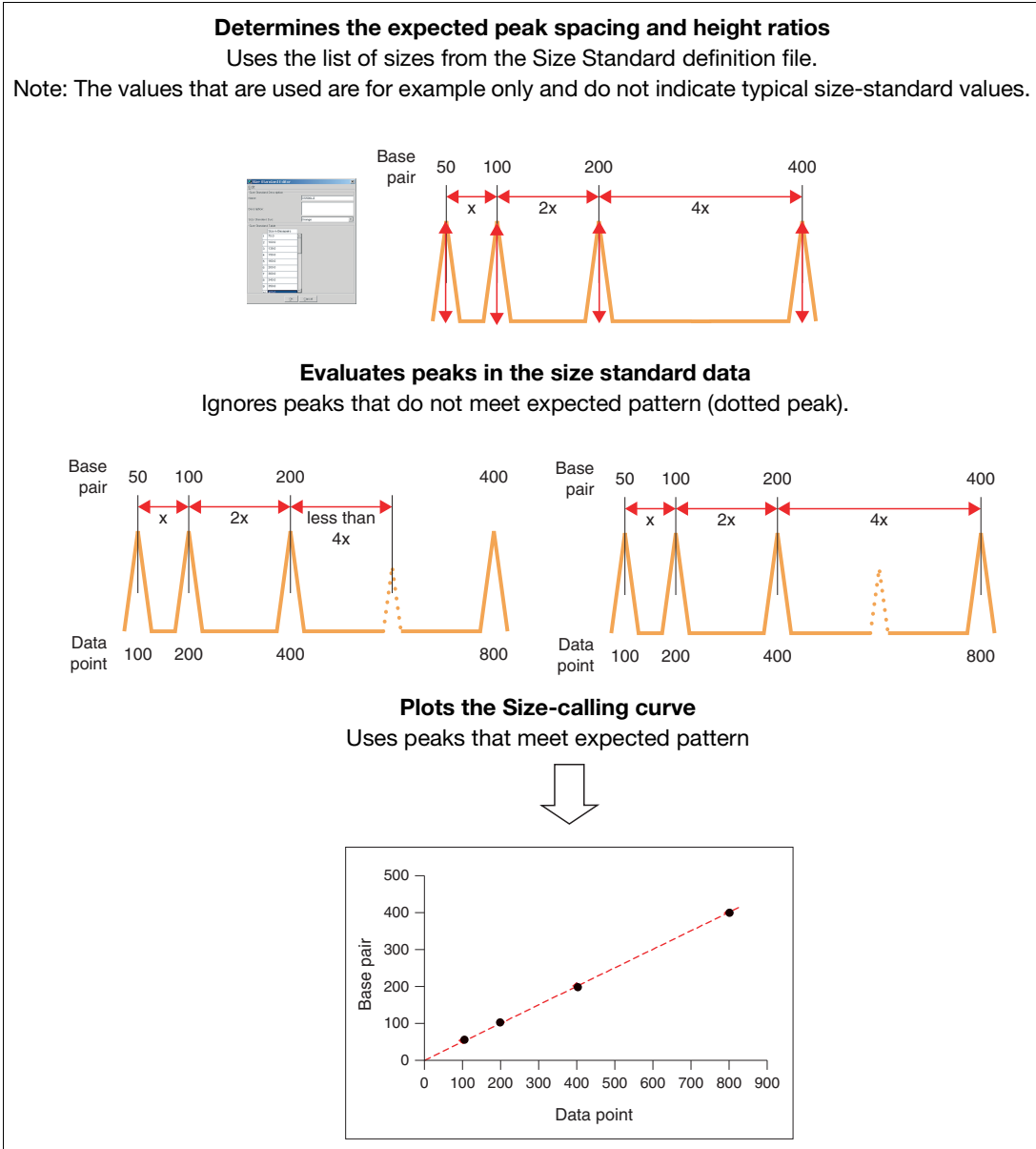


Figure 11 Size matching and size-calling curve generation

## Size-Calling Curve Generation and Size Calling

### Sizing:

- Size matching
- **Size-calling curve generation**
- **Size calling**
- Sizing quality

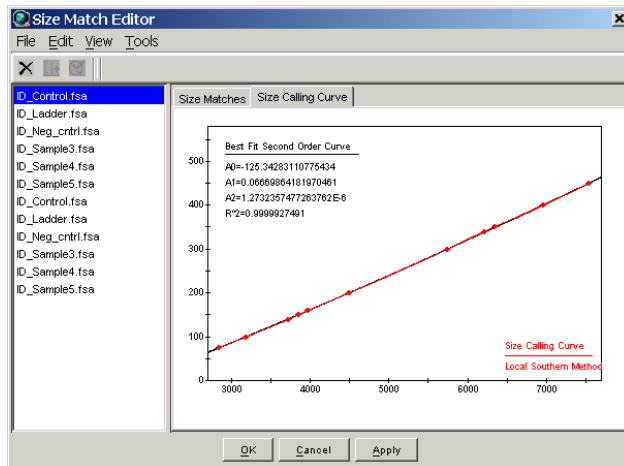
To generate the size-calling curve, the software plots the actual data points of the size standard against the expected size of each size standard peak. The size-calling method determines how the size-calling curve is generated and used to size each sample.

During size matching and size calling:

- Two size-calling curves are generated for each sample:
  - Black: A best-fit second order curve, regardless of the size-calling method that is selected.
  - Red: A curve based on the size-calling method that is selected in the analysis method.
- The data points of non-size-standard peaks are plotted against the size-calling curve
- Peaks are sized according to the size-calling method that is selected in the analysis method

Size-calling methods are:

- Local Southern
- Global Southern
- Least Squares (2nd Order or 3rd Order)
- Cubic Spline Interpolation



## Size Calling – Local Southern Method

- Sizing:
- Size matching
  - Size-calling curve generation
  - **Size calling**
  - Sizing quality assessment

The Local Southern method (Figure 12) determines the sizes of fragments by using the reciprocal relationship between fragment length and mobility, as described by E. M. Southern (1979).

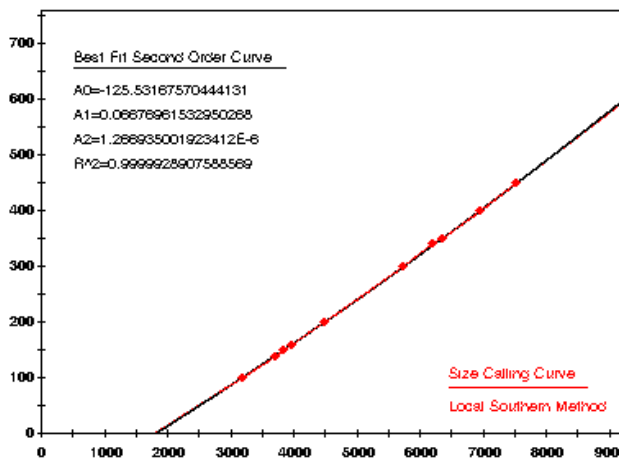


Figure 12 Local Southern method

**IMPORTANT!** Do not use the Local Southern size-calling method for AmpF $\mathcal{L}$ STR<sup>®</sup> MiniFiler<sup>™</sup> PCR Amplification Kit data. The Local Southern method requires two data points below the smallest fragment for accurate sizing. Because of the small size of amplicons generated using the MiniFiler<sup>™</sup> kit, the GeneScan<sup>™</sup> 500 LIZ<sup>®</sup> Size Standard 50 bp peak may be hidden by the primer peak, and allele 5 in the D16S539 marker can size at <75 bp. Use the 3rd Order Least Squares size-calling method described on [page 22](#).

### Local Southern Method Equation

The equation attempts to describe the reciprocal relationship between the mobility,  $m$ , and the length,  $L_0$ , of the standard fragments.

$$L = [c/(m - m_0)] + L_0$$

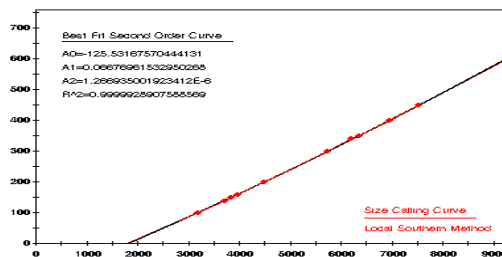
## How The Local Southern Method Works

This method, which is similar to the Cubic Spline method (page 24), uses the four fragments closest in size to the unknown fragment to determine a best-fit line value. Only the region of the size standard near the fragment of unknown length is analyzed.

**Note:** Size estimates may be inaccurate if any of the standard fragments run anomalously.

In the Local Southern method (Figure 13):

1. The fitting constants of the curve are calculated for each group of three neighboring points on the standard.
2. A curve is then created using three standard points (two points below and one point above the fragment), then a fragment size is determined.
3. Another curve is created using an additional set of three points (one point below and two points above the fragment), to assign another value.
4. The two curves are averaged to determine the unknown fragment length.



To size call fragments between b and c:

- Fit curve through a, b, and c
- Fit curve through b, c, and d
- Average curves between b and c

Figure 13 Local Southern method

## Size Calling – Global Southern Method

- Sizing:
- Size matching
  - Size-calling curve generation
  - **Size calling**
  - Sizing quality assessment

The Global Southern method (Figure 14) is similar to the Least Squares method (page 22) 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 on that line to calculate the fragment values.

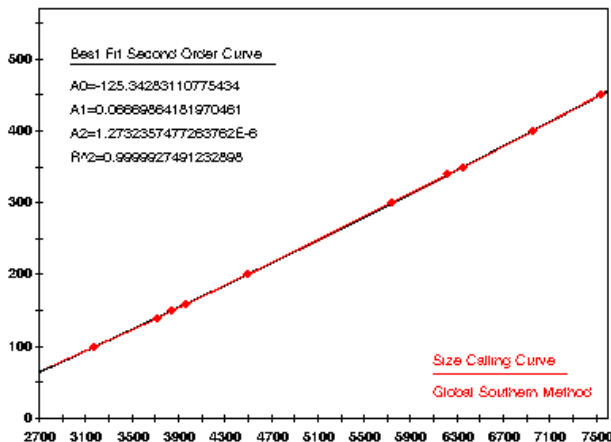


Figure 14 Global Southern method

### Global Southern Method Equations

Equation	Description
$L = [c / (m - m_0)] + L_0$	Attempts to describe the reciprocal relationship between the mobility, m, and the length, L <sub>0</sub> , of the standard fragments.
$\sum i \{ L_i - [c / ((m_i - m_0) + L_0)] \}^2$	The fitting constants L <sub>0</sub> , m <sub>0</sub> , and c are calculated by a least-squares fit to minimize the left side quantity.

### How the Global Southern 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. For best results, use a standard that brackets all the fragments of interest.

## Size Calling – Least Squares Method

### Sizing:

- Size matching
- Size-calling curve generation
- **Size calling**
- Sizing quality assessment

Both Least Squares methods (2nd-Order and 3rd-Order) use regression analysis to build a best-fit size-calling curve. This curve:

- Produces the minimum additive distance from the curve to the plotted data points.
- Compensates for any fragments that may run anomalously.

Consequently, this method typically results in the least amount of deviation 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 data points to produce a sizing curve based on Multiple Linear Regression.

### Using 3rd Order for MiniFiler™ Kit Data

A 3rd-order curve (which uses a higher polynomial degree and captures more of the peak structure) provides the most flexibility when generating best-fit curves for sizing samples with anomalously migrating fragments.

### Examples

[Figure 15 on page 23](#) and [Figure 16 on page 24](#) show 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.

To generate the Least Squares curve, the software:

- Plots the known fragment sizes (bp) versus data points.
- Generates a best-fit curve using regression analysis.



- Applies the following calculation to determine the size in data points of the unknown fragments:

$$Y = Ax^3 + Bx^2 + Cx + D$$

where:

Y = size (bp)

x = datapoint

A = First order coefficient

B = Second order coefficient

C = Third order coefficient

D = Zeroth Coefficient or constant

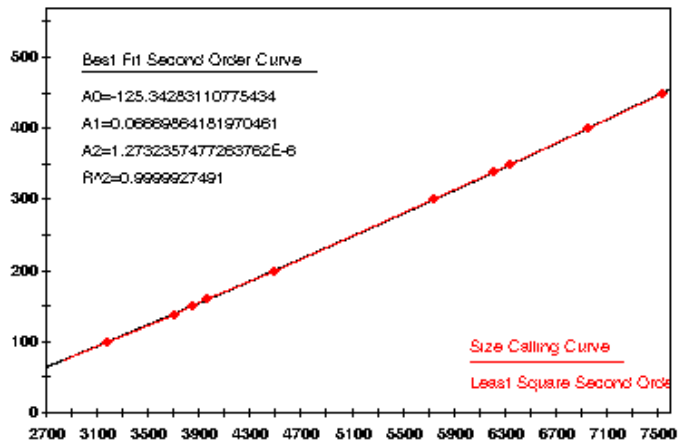


Figure 15 2nd-Order Least Squares size-calling curve (quadratic)

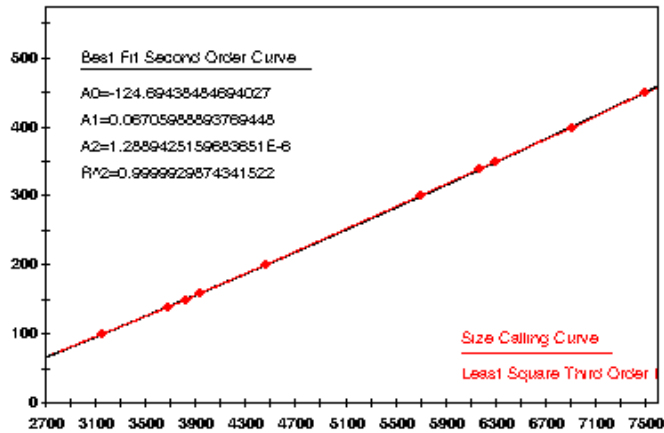


Figure 16 3rd-Order Least Squares size-calling curve (cubic)

## Size Calling – Cubic Spline Interpolation Method

### Sizing:

- Size matching
- Size-calling curve generation
- **Size calling**
- Sizing quality assessment

The Cubic Spline method (Figure 17), which is similar to the Local Southern method (page 19), forces the sizing curve through all the known points of the selected size standard. Although this enforcement produces exact results for the values of the standards themselves, it does not compensate for standard fragments that may run anomalously.

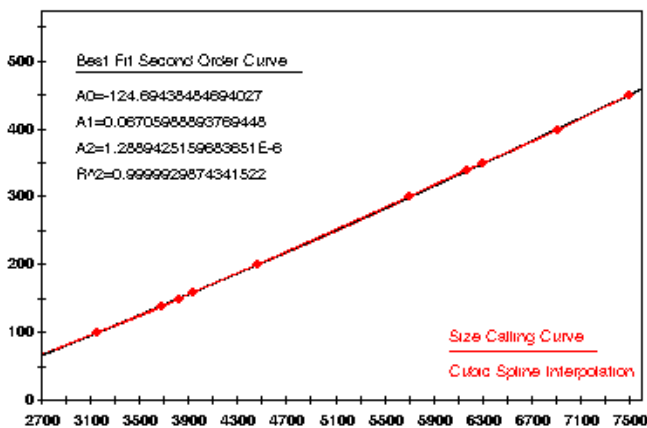


Figure 17 Cubic Spline Interpolation method

## 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 degree of sizing accuracy error.


---

## Sizing Quality Assessment

Sizing:

- Size matching
- Size-calling curve generation
- Size calling
- **Sizing quality assessment**

The last step in sizing is the sizing quality assessment. Sizing quality is discussed in [“Sizing Quality Assessment” on page 43](#).

The software does not genotype samples with  sizing quality.

## Allelic Ladder Genotyping

### Allelic ladder genotyping:

- Allele matching
- Allelic ladder quality assessment
- Bin offsetting

After sizing quality is determined, the following occur:

- Allele matching
- Allelic ladder quality assessment
- Bin offsetting

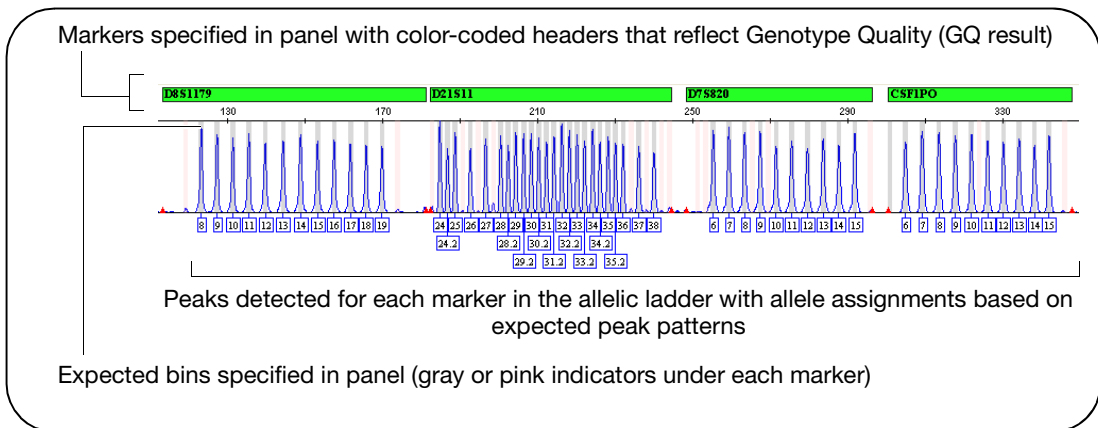
## Allele Matching

### Allelic ladder genotyping:

- **Allele matching**
- Allelic ladder quality assessment
- Bin offsetting

During allele matching, the software assigns allele labels to allelic ladder peaks by matching peaks with the bins that are specified in the panel. The matching algorithm uses ratio-matching, based on relative distance and height of neighboring peaks.

Figure 18 is an example of the blue dye plot in the AmpF $\Phi$ STR<sup>®</sup> Identifiler<sup>®</sup> PCR Amplification Kit allelic ladder.



**Figure 18** Identifiler<sup>®</sup> kit allelic ladder blue dye

## Allelic Ladder Quality Assessment

Allelic ladder genotyping:

- Allele matching
- **Allelic ladder quality assessment**
- Bin offsetting

After allele matching, allelic ladders are evaluated to determine whether they are used to create bin offsets, described on [page 27](#).

Allelic ladder quality assessment is discussed in “[Allelic Ladder Quality Assessment](#)” on [page 45](#).

## Bin Offsetting

Allelic ladder genotyping:

- Allele matching
- Allelic ladder quality assessment
- **Bin offsetting**

Bin offsetting evaluates the alleles found in allelic ladder samples and adjusts bins before allele-calling. Marker ranges are also adjusted based on the bin offsets.

### AmpF $\lambda$ STR<sup>®</sup> Allelic Ladder and Bin Sets

The allelic ladder for each AmpF $\lambda$ STR<sup>®</sup> kit contains the most common alleles for each locus and is used to generate the reference sizes in the bin sets provided with the software.

Reference sizes consider:

- The actual size (bp) of DNA fragments.
- 3' A nucleotide addition.
- Mobility shifts related to the non-nucleotide linkers in the allelic ladder. Non-nucleotide linkers enable reproducible positioning of the alleles to facilitate inter-locus spacing. (Grossman *et al.*, 1994 and Baron *et al.*, 1996)

## Physical and Virtual Bins

The bin set provides reference allele sizes for:

- Alleles physically present in the allelic ladder (physical bins).
- Alleles that are not present in the allelic ladder (virtual bins), but have either been reported in STRbase ([www.cstl.nist.gov/div831/strbase](http://www.cstl.nist.gov/div831/strbase)) or discovered during developmental validation of a particular chemistry at Applied Biosystems, and can be genotyped by reference to the alleles present in the ladder.

Figure 19 illustrates a bin set with eight physical bins and six virtual bins.

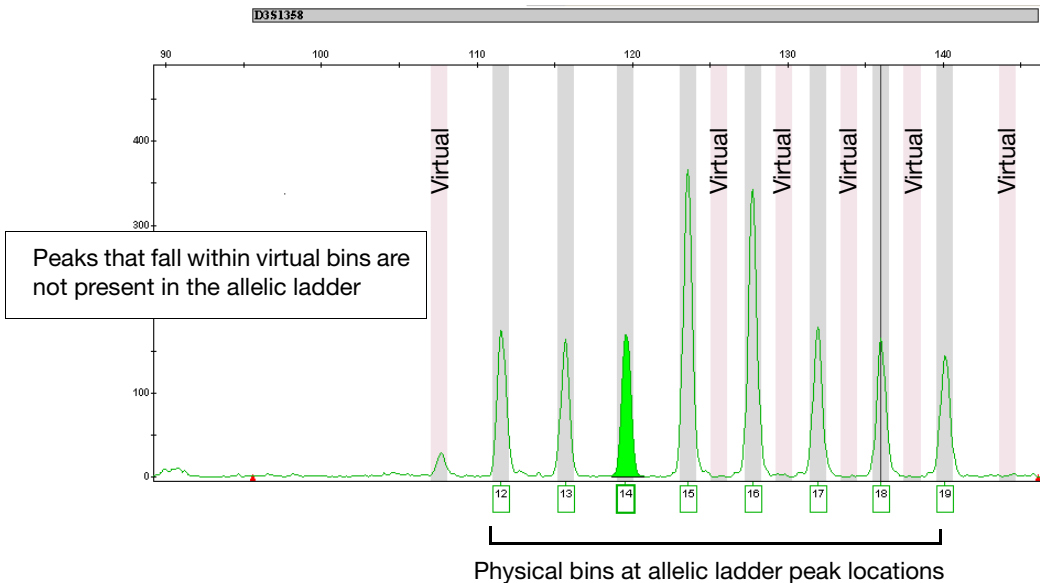




Figure 19 Bin set with physical and virtual bins

**Note:** The GeneMapper® *ID-X* Software allows you to display virtual bins in pink. For information, see the GeneMapper® *ID-X* Software Help.

## How the Software Calculates and Applies Bin Offsets

When analyzing allelic ladder samples, the GeneMapper® *ID-X* Software offsets the reference sizes defined in the bin set relative to the actual sizes obtained for the alleles in the allelic ladder.

To determine offsets, the software:

1. Performs an allelic ladder quality assessment (described on [page 45](#)).
2. Uses  CGQ allelic ladders with  SQ only. CGQ is described on [page 54](#). SQ is described on [page 43](#).
3. Identifies the size of a physical allele in the allelic ladder. Virtual alleles are not used to calculate offsets.  
If more than one allelic ladder sample is present in the run folder, determines the offset for each allele in each of the allelic ladder samples, then averages the offsets for the allele.
4. Offsets the corresponding bins for the physical alleles by the difference between the observed size for the allele in the run data and the reference size for the allele.
5. Offsets the virtual bins. To offset the virtual bins, the software uses the offset from the closest physical bin or virtual bin to the left of the bin in question. If the first bin is a virtual bin, the software uses the offset from the first physical bin (counting from left to right).

**6.** Offsets the:

- Marker Range Min Size value by the calculated bin offset of the first physical bin of the allelic ladder in the marker.
- Marker Range Max Size value by the calculated bin offset of the last physical bin of the allelic ladder in the marker.

For more information, see [“Displaying Offset Marker Margins” on page 33](#).

---

**IMPORTANT!** The software displays the *reference* marker range in two places: in the Samples plot status bar when you click the marker header above the plot and in the Panel Manager. To display the *offset* marker ranges, examine the value that is displayed in the Samples plot status bar when you click the red marker indicators on the x-axis of the Samples plot.

---

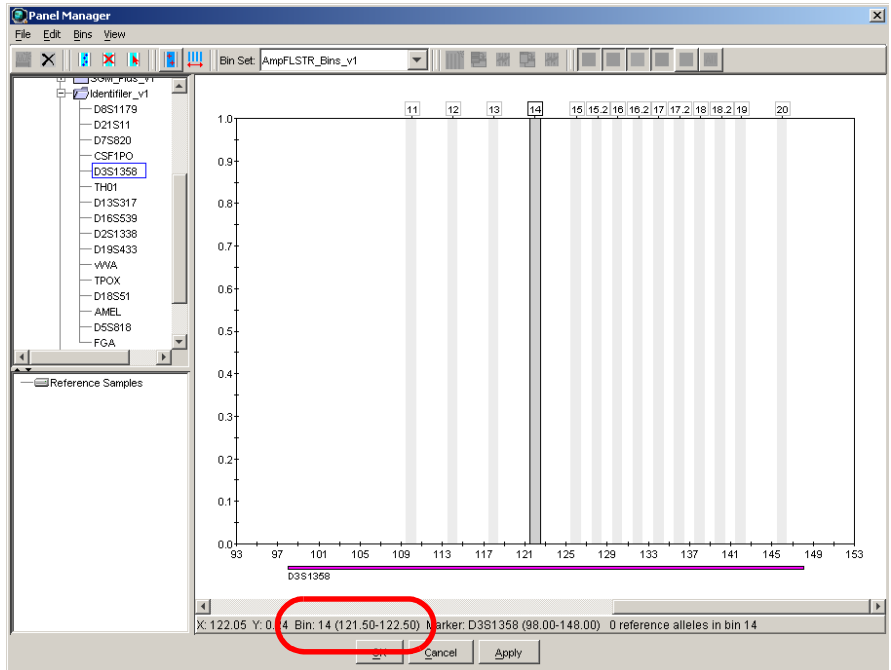
**7.** Repeats the process for each physical allele in all samples with a Sample Type of Allelic Ladder.

**Example: One Allelic Ladder Is Present In a Run Folder**

A reference size in the bin set for an allele is 122 bp. The observed size is 119.56 bp. The offset value is calculated as  $119.56 - 122 = -2.44$ . The bin for this allele is shifted down 2.44 base pairs to 119.56 ( $\pm 0.5$  bp).

[Figure 20 on page 31](#) shows the reference size (bin) for allele 14 in the D3S1358 marker before analysis.

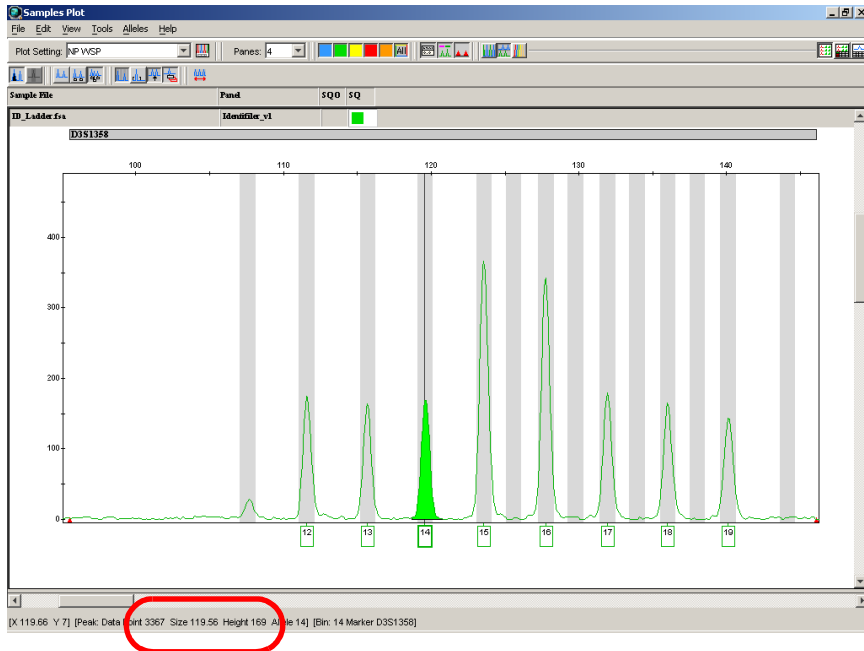




Reference allele size (position of the bin)  $122 \pm 0.5$  bp

Figure 20 Bin assignment before offsetting

Figure 21 on page 32 shows that after analysis of the allelic ladder sample, the bin is offset based on the observed size of the allele in the allelic ladder.



Offset bin  $119 \pm 0.5$  bp

**Figure 21** Bin assignment after offsetting

**Example: More Than One Allelic Ladder Sample Is Present In a Run Folder**

A reference size in the bin set for an allele is 122 bp. The observed size is 119.56 bp for one allelic ladder sample and 119.50 bp for another allelic ladder sample. The offset for the each allelic ladder sample is calculated  $(119.56 - 122 = -2.44)$  and  $(119.50 - 122 = -2.50)$ . The software averages the two offsets  $[(-2.44 + -2.50)/2 = -2.47]$ . The bin for this allele is shifted down 2.47 base pairs to  $119.53 (\pm 0.5$  bp).

**Example: How Offsets Are Applied To Virtual Alleles**

Virtual allele 11 in [Figure 21](#) above uses the same offset value as allele 12. The sizes between these two alleles differ by 4 bp. Specifying a size for allele 11 that is 4 bp shorter than allele 12 is a reasonable estimate because alleles 11 and 12 differ by a single repeat unit (4 bp).

## Displaying Offset Marker Margins

When bins are offset, marker margins are also offset. To display the shifted marker margins, click the red marker indicators in the Samples plot (Figure 22).

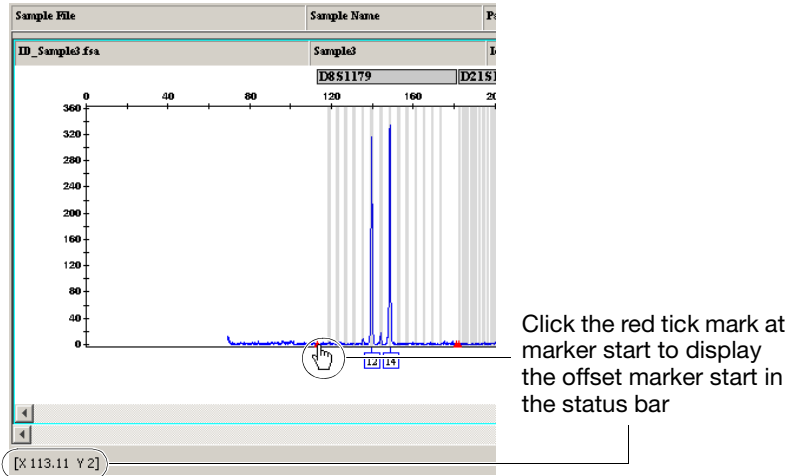


Figure 22 Displaying offset marker margins

**Note:** Reference marker margins that are defined in the panel are displayed when you click the marker header in the Samples plot and in the Panel Manager (Figure 23).

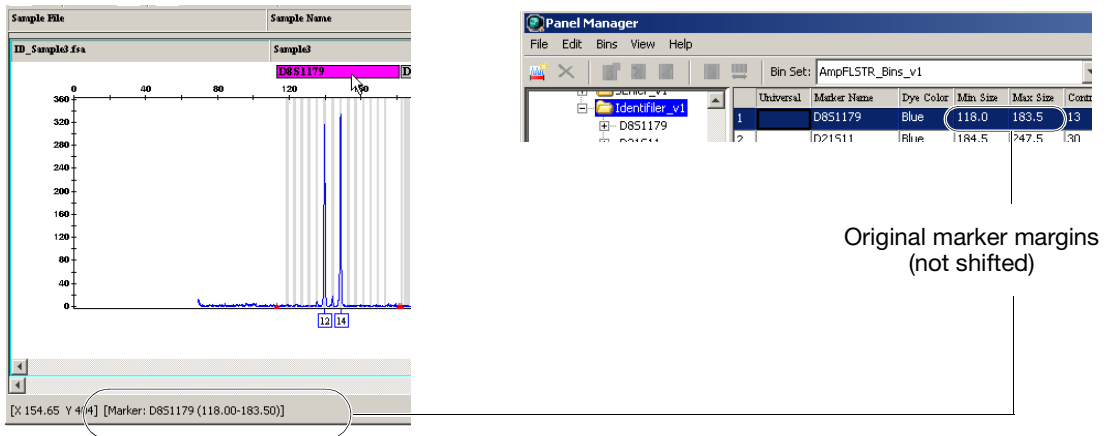


Figure 23 Original marker margins are displayed when you click the marker header in the Samples plot and in the Panel Manager

## Sample Genotyping

### Sample genotyping:

- Allele calling and filtering
- Genotype and sample and quality assessment

After allelic ladders are detected and sized, sizing quality is determined, then allelic ladders are genotyped.

After allelic ladders are genotyped, samples are detected and sized, sizing quality is determined, then samples are genotyped.

During genotyping of samples, the GeneMapper® *ID-X* Software:

- Performs allele-calling and filtering.
- Assesses genotype quality.
- Assesses sample quality.

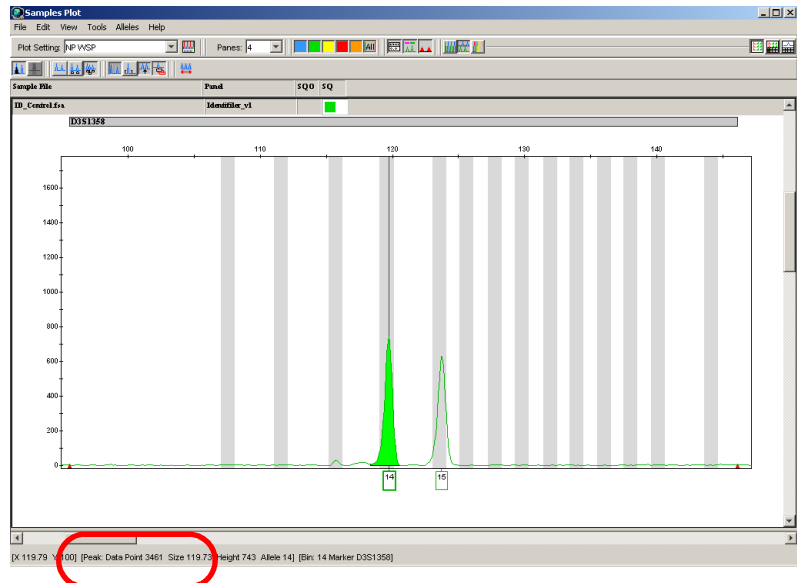
## Allele Calling and Filtering of Samples

### Sample genotyping:

- **Allele calling and filtering**
- Genotype and sample and quality assessment

During allele calling of samples, the software:

- Labels allele peaks in unknown samples:
  - Compares sizes of the unknown peaks to the offset bins.
  - Assigns the relevant allele label if a peak falls within one of the defined bins. [Figure 24 on page 35](#) is an example of a 14 allele in D3S1358.
  - Assigns an OL (Off-Ladder) label if a peak falls outside of the defined bins.
- Applies filters to eliminate peaks from consideration as allele peaks:
  - Allele Tab analysis method settings (Global Cut-off value, Minus A, Global Stutter, Amelogenin Cutoff, and Range Filter)
  - Marker-specific stutter ratios if specified in the Analysis Method (Allele tab) and defined in the panel



Reference allele size (position of the bin)  $119 \pm 0.5$  bp

**Figure 24 Allele calling**

Notice the size of this allele is 119.73, which is within the shifted bin calculated previously ( $119.53 \pm 0.5$  bp, see [Figure 21 on page 32](#)). Therefore, this peak is labeled as a 14.

**Off Ladder (OL)  
Allele Calls**

If a peak falls outside one of the defined bins, the software labels it OL (Off Ladder).

OL calls may be caused by:

- The presence of a microvariant allele at a particular locus
- Sample migration anomalies and artifacts

**Allele Naming  
Convention**

Allele names are assigned based on the number of complete four-base-pair repeat-units and partial repeat-units:

- Complete four base pair repeat units are designated by an integer.
- Variant alleles that contain a partial repeat are designated by a decimal followed by the number of bases in the partial repeat. For example:


- An FGA 26.2 allele contains 26 complete repeat units and a partial repeat unit of two base pairs.
- A D21S11 allele 33.1 contains 33 complete repeat units and one nonconsensus base pair.
- A D21S11 29.3 allele contains 29 complete repeat units and a partial 3-bp unit (Moller et al., 1995; Gill et al., 1997).

## Genotype and Sample Quality Assessment

### Sample genotyping:

- Allele calling and filtering
- **Genotype and sample and quality assessment**

To determine sample quality, the software:

- Applies user-defined and software-defined thresholds to generate marker-level quality value flags (described in [“Marker-Level Quality Assessment” on page 47](#)) and sample-level quality value flags (described in [“Sample-Level Quality Assessments” on page 53](#)).
- Applies GQ weighting from  marker-level quality value flags to determine the Genotype quality (GQ) (described in [“Sample-Level Quality Assessments” on page 53](#)).
- Uses GQ to determine the sample-level CGQ composite genotyping quality assessment (described in [“CGQ \(samples\)” on page 54](#)).

## 2

# GeneMapper® *ID-X* Software Quality Value System

This chapter covers:




- Overview of the Quality Value System . . . . .38
- Overview of the Quality Value System Checks and Assessments. . . . .40
- Analysis Requirement Checks. . . . .41
- Sizing Quality Assessment . . . . .43
- Allelic Ladder Quality Assessment. . . . .45
- Marker-Level Quality Assessment . . . . .47
- Genotype Quality Assessment. . . . .51
- Sample-Level Quality Assessments . . . . .53

## Overview of the Quality Value System

The GeneMapper® *ID-X* Software provides a quality value system and a set of streamlined data review tools and features that can be optimized and validated for use in both expert systems and traditional data analysis workflows.

The GeneMapper® *ID-X* Software quality value system:

- Assesses the quality of allelic ladders before analysis and does not consider low-quality allelic ladders for genotyping.
- Assesses the quality of data at the sample and marker level using process quality values (PQVs).
- Can be used in an optimized and validated expert system or traditional data analysis workflow to quickly identify data quality issues and aid in interpretation of samples that do not meet all thresholds.
- Can be used in an optimized and validated expert system workflow to quickly segregate samples that require manual review from those that do not.

The PQV results of the quality assessment are displayed as color-coded flags:   . The color of the flag depends on software-specified thresholds or user-defined thresholds set in the analysis method.

### Optimizing and Validating an Expert System

Before using any software as an expert system, optimize and validate the thresholds for each AmpF $\Phi$ STR kit and instrument platform combination by processing a variety of samples that challenge each of the quality flags.

---

**IMPORTANT!** Different kit/instrument combinations may require different thresholds.

---



## Displaying Numeric SQ and GQ Values

By default, the GeneMapper *ID-X* Software displays Pass ■, Check-▲, or Low Quality ● for Sizing Quality (SQ) and Genotyping Quality (GQ) to represent the numeric score of the assessment.

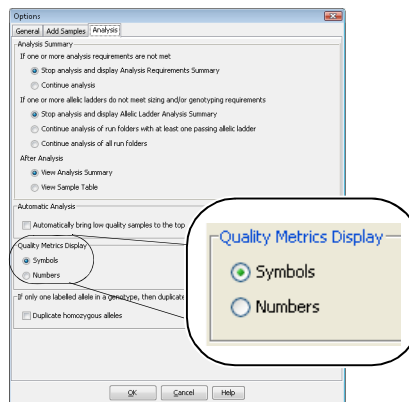
For example, the Sizing Quality (SQ) yields a value between 1 and 0. Based on the SQ threshold settings of the SQ & GQ Settings tab, the software displays Pass ■, Check-▲, or Low Quality ● to indicate the SQ result.

You can alternatively display numeric values for SQ and GQ.

Sizing Quality Representation	Samples Table Example																		
Symbols (default)	<table border="1"> <thead> <tr> <th>Run Name</th> <th>SFNF</th> <th>SNF</th> <th>OS</th> <th>SQ</th> <th>WELLQ</th> </tr> </thead> <tbody> <tr> <td>DGB_SNPlex_</td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> </tr> <tr> <td>DGB_SNPlex_</td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> </tr> </tbody> </table>	Run Name	SFNF	SNF	OS	SQ	WELLQ	DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>
Run Name	SFNF	SNF	OS	SQ	WELLQ														
DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>														
DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>														
Numbers	<table border="1"> <thead> <tr> <th>Run Name</th> <th>SFNF</th> <th>SNF</th> <th>OS</th> <th>SQ</th> <th>WELLQ</th> </tr> </thead> <tbody> <tr> <td>DGB_SNPlex_</td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td>1.0</td> <td>1.0</td> </tr> <tr> <td>DGB_SNPlex_</td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td><span style="color: green;">■</span></td> <td>1.0</td> <td>1.0</td> </tr> </tbody> </table>	Run Name	SFNF	SNF	OS	SQ	WELLQ	DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	1.0	1.0	DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	1.0	1.0
Run Name	SFNF	SNF	OS	SQ	WELLQ														
DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	1.0	1.0														
DGB_SNPlex_	<span style="color: green;">■</span>	<span style="color: green;">■</span>	<span style="color: green;">■</span>	1.0	1.0														

To display numeric values for SQ and GQ:

1. Select **File ▶ Project Options**, then select the **Analysis** tab.
2. In the Quality Metrics Display settings, select **Numbers**.



3. Click **OK** to apply the settings.

## Overview of the Quality Value System Checks and Assessments

The GeneMapper *ID-X* Software quality value system performs the following checks and assessments:

- **Analysis requirements checks:**

Before analysis starts, identifies any conditions that may prevent analysis or cause unexpected results. If one or more analysis requirements are not met, the ARNM (Analysis Requirements Not Met) PQV is set to ▲. If Project Options are set accordingly, the Analysis Requirements Summary is also displayed (described in GeneMapper® *ID-X* Software Help).

- **Allelic ladder genotyping:**

- **Sizing quality assessment** – Evaluates the quality of the size standard profile in each allelic ladder.

- **Allelic ladder quality assessment** – Evaluates allelic ladder quality. Also determines if an allelic ladder is used for creating bin offsets. If the software finds one or more allelic ladders that do not meet sizing and genotyping requirements and Project Options are set accordingly, the Allelic Ladder Analysis Summary is displayed (described in GeneMapper® *ID-X* Software Help).

- **Marker-level quality assessment** – Evaluates labeled peaks within each marker. The SPK and OS PQVs contribute to the allelic ladder GQ assessment if spikes and/or off-scale fluorescence signals are present in the allelic ladder (and Spike Detection is Enabled for the Allelic Ladder and the Allelic Ladder GQ Weighting for Spikes and Off-scale is set to 1[on]. If Spike Detection is Disabled for the Allelic Ladder, spikes will not be detected or labeled.) Other PQVs do not affect allelic ladder GQ.


- **Sample-level quality assessment** – Evaluates the quality of the entire sample.

- **Sample genotyping:**







- **Sizing quality assessment** – Evaluates the quality of the size standard profile in each sample.

- **Marker-level quality assessment** – Evaluates labeled peaks within each marker. Contributes to the genotype quality assessment.
- **Sample-level quality assessment** – Evaluates the quality of the entire sample.
- **Genotype quality assessment** – Evaluates the quality of each marker in a sample. Contributes to the overall composite genotype quality assessment.

## Analysis Requirement Checks



The analysis requirements checks are performed before data analysis starts. If one or more analysis requirement checks is , analysis may not occur or may not give the expected results (for example, analyzing without a size standard or allelic ladder affects sizing quality or genotyping quality).

Analysis requirements checks that are displayed in the Samples table before the analysis starts




Acronym	Full Name	Function/Rule(s)	Flag	
SFNF	Sample File Not Found	Indicates if the software finds the sample (.fsa or .hid) files in the original location from which they were added to the project.  Select <b>Edit ▶ Define New Sample File Path</b> to re-associate the samples file with the project.		Sample file found
				Sample file not found
MNF	Matrix Not Found	Indicates if the software finds in the database the matrix file that is associated with the sample.		Matrix file found
				Matrix file not found
SNF	Size Standard Not Found	Indicates if the software finds in the database the size standard definition associated with the sample.		Size standard definition found
				Size standard definition not found

**Note:** The SFNF, MNF, and SNF flags are included in the ARNM flag described on [page 42](#). To conserve space, display only the ARNM column in the Samples table.




Analysis requirements checks that are displayed in the Analysis Requirements Summary after the analysis starts

Acronym	Full Name	Function/Rule(s)	Flag Indicator	
ARNM	Analysis Requirement Not Met	<p>Indicates if all analysis requirements are met. These requirement checks are performed when analysis is started, before actual analysis begins:</p> <ul style="list-style-type: none"> <li>• Sample File Not Found</li> <li>• Analysis Method Not Selected</li> <li>• Analysis Method Not Found In The Database</li> <li>• Panel Not Selected</li> <li>• Panel Not Found in the Database</li> <li>• Binset Not Selected</li> <li>• Binset In Analysis Method does not match Binset selected in the Panel Manager</li> <li>• Size Standard Not Selected</li> <li>• Size Standard Not Found In Database</li> <li>• Size Standard Dye Color Is Not Present In The Sample Dye Set</li> <li>• Matrix Not Selected</li> <li>• Matrix Not Found or Contains Invalid Data</li> <li>• No Allelic Ladder Selected In Run Folder</li> <li>• GMID V3.x Analysis Method Selected</li> <li>• Basic Or Classic Size Standard Selected</li> <li>• SNP Panel Selected</li> </ul> <p><b>Note:</b> The ARNM PQV includes the SFNF, MNF, and SNF PQVs described on the previous page. To conserve space, display only the ARNM column in the Samples table.</p>		All analysis requirements are met.
				One or more analysis requirements are not met.



## Sizing Quality Assessment

The quality value system evaluates the quality of the size standard profile within each sample (SQ) and allows you to flag size standards with poor peak resolution. You can adjust the ranges that correspond to the Pass , Check , and Low Quality  ranges in the analysis method (SQ & GQ Settings).

Sizing Quality that is displayed in the Samples Table after the analysis is complete

Acronym	Full Name	Function/Rule(s)	Flag Indicator	
SQ	Sizing Quality	<p>Evaluates the similarity between the fragment pattern for the size standard dye that is specified in the size standard definition and the actual distribution of size standard peaks in the sample, calculates an interim SQ (a value between 0 and 1) as described in <a href="#">“Sizing” on page 16</a>, then applies the broad peak weighting that is specified in the analysis method.</p> <p><b>Weighting</b></p> <p>The Broad Peak (BD) threshold and SQ Broad Peak weighting that is specified in the analysis method affect the SQ. To determine the final SQ value, the software:</p> <ul style="list-style-type: none"> <li>Evaluates peak widths in the sample in the dye color specified in the size standard definition.</li> <li>If the width of any peak in the sizing range exceeds the broad peak threshold, applies the SQ weighting: Interim SQ × (1 – SQ weighting)</li> </ul>		<p>SQ is within the user-defined Passing range (Default = 0.75 to 1.0)</p>
				<p>SQ is between the user-defined Passing range and Low Quality range (Default = 0.26 to 0.74)</p>
				<p>SQ is within the user-defined Low Quality range (Default = 0.0 to 0.25)</p>

### Sizing Quality that is displayed in the Samples Table after the analysis is complete

Acronym	Full Name	Function/Rule(s)	Flag Indicator
SQ (continued)	Sizing Quality	<p><b>Example</b></p> <p>A size standard peak exceeds BD threshold, has an Interim SQ = 0.8, and SQ weighting = 0.5. The final SQ is:</p> $SQ = 0.8 \times (1 - 0.5) = 0.4$ <p>Based on the default SQ range, this SQ is .</p> <p><b>Note:</b> Broad Peak (BD) threshold and SQ Broad Peak weighting are user-defined.</p> <p><b>Note:</b> The GeneMapper® ID-X Software does not genotype samples with  SQ.</p>	

## Allelic Ladder Quality Assessment

The quality value system performs an allelic ladder quality assessment to determine if a ladder is used in genotyping (to create bin offsets, described on [page 27](#)).


Allelic ladder quality is reported per marker by the GQ (Genotyping Quality is described on [page 51](#)) and per sample by the CGQ (Genotyping Quality is described on [page 54](#)).

### Quality Rules


Allelic ladder samples are analyzed before all other samples. An allelic ladder sample must have a   SQ and a   CGQ to be used for creating bin offsets. For an allelic ladder to have a   CGQ, all the markers within the allelic ladder must pass the following rules:

Rule	Description
1	All ladder alleles specified in the panel used to analyze are detected.
2	In each marker, the peak height ratio of the first and second peak is greater than 50%. This rule eliminates allelic ladders if the stutter peak before the first true allele peak is labeled as an allele.
3	No spikes are detected above 20% (default) of the highest allele peak in the same dye color within the extended marker range. <b>Note:</b> When spike detection is enabled for Allelic Ladders, spike detection is performed within each extended marker range (no gaps are present between markers; the end point of each marker is extended past the marker definition in the panel to the beginning of the next marker). <b>Note:</b> When spike detection for allelic ladders is set to Disable, spikes will not be detected or labeled. <b>Note:</b> The Allelic Ladder Spike Cut-off value is user-definable in the Peak Quality tab of the analysis method.
4	The peak height ratio between the lowest and highest peak is equal to or greater than 15%.
5	In each marker, the base pair spacing between any two ladder alleles specified in the panel used to analyze is within the expected range.
6	No off-scale (OS) fluorescent signal is detected within each extended marker range. <b>Note:</b> The Allelic Ladder GQ Weighting for Off-scale is user-definable.

**Effect on GQ** Allelic ladder quality rules are evaluated at the marker level and affect GQ results.

If any of the rules do not pass, the GQ (described on [page 51](#)) for the allelic ladder marker is set to .

---

**Note:** If rule 3 does not pass and the Allelic Ladder GQ Weighting for Spikes is set to 0 (off in the SQ & GQ Settings tab in the analysis method), the GQ for the allelic ladder marker is set to , even if spikes are present.

---



Other than SPK, no other individual marker-level PQV flags (described on [page 47](#)) affect the GQ of allelic ladder samples. Therefore, they do not affect the allelic ladders used for genotyping.







## Marker-Level Quality Assessment

Marker-level quality assessment indicates the quality of each marker in a sample.







Marker-level quality assessment evaluates only labeled peaks within a defined marker size range defined in the panel and can be weighted to contribute to the calculation of the GQ (Genotype Quality).

**Note:** Gray PQVs ( ) indicate that at least one label in the marker was edited or the GQ was overridden.







Marker-level quality flags that are displayed in the Genotypes table after the analysis is complete

Acronym	Full Name	Function/Rule(s)	Flag Indicator	
OS	Off-scale	Indicates if any fluorescence signal exceeds the detection threshold of the instrument.  <b>Note:</b> The detection limit varies on different instruments. The OS quality flag indicates if the CCD camera is saturated at any data point within the analysis range. It is not based on an absolute RFU value (peak height).		No fluorescence signal exceeds the detection limit of the instrument.
				One or more fluorescence signals exceed the detection limit of the instrument.
BIN	Out of Bin Allele	Indicates if labeled peaks do not fall inside bins. These peaks are labeled with OL (off ladder). For information of bin offsetting, see <a href="#">page 27</a> .		No labeled peaks are detected outside an offset bin.
				One or more labeled peaks are detected outside an offset bin.





Marker-level quality flags that are displayed in the Genotypes table after the analysis is complete (*continued*)

Acronym	Full Name	Function/Rule(s)	Flag Indicator	
PHR	Peak Height Ratio	Indicates if the peak height ratio between the lowest and highest peak is less than the Min Peak Height Ratio defined in the analysis method.		The peak height ratio calculated between the lowest and highest peaks is equal to or greater than the Min Peak Height Ratio threshold.
				The peak height ratio calculated between the lowest and highest peaks is less than the Min Peak Height Ratio threshold.
			N/A	Only one peak is labeled.
MPH	Max Peak Height	Indicates if any peak heights (RFU) exceed the Max Peak Height (RFU) value set in the analysis method.		No peaks exceed the Max Peak Height (RFU) value (all peak heights are less than or equal to the threshold).
				One or more peaks exceed the Max Peak Height (RFU) value.
			N/A	No peaks are detected or all peaks are off-scale (OS).
LPH	Low Peak Height	Indicates if any peak heights (RFU) are below the following thresholds set in the analysis method: <ul style="list-style-type: none"> <li>• Homozygous Min Peak Height</li> <li>• Heterozygous Min Peak Height</li> </ul>		All detected peak heights are equal to or greater than the thresholds.
				One or more peak heights are below the thresholds.

Marker-level quality flags that are displayed in the Genotypes table after the analysis is complete (*continued*)

Acronym	Full Name	Function/Rule(s)	Flag Indicator	
AN	Allele Number	Indicates if the software detects: <ul style="list-style-type: none"> <li>No alleles <i>or</i></li> <li>More than the Max Expected Alleles set in the analysis method (Peak Quality tab) <i>or</i></li> <li>No X allele in Amelogenin</li> </ul>		At least one allele and not more than the Max Expected Alleles are detected or X allele for Amelogenin is detected.
				No alleles, more than the Max Expected Alleles, or no X allele for Amelogenin is detected.
BD	Broad Peak	Indicates if the width of any peak exceeds the Max Peak Width (half height in basepairs) defined in the analysis method (Peak Quality tab).  Broad peaks may be a result of dye artifacts or poor resolution. Peaks with poor resolution: <ul style="list-style-type: none"> <li>May be caused by electrophoresis issues</li> <li>May not be true DNA peaks</li> </ul>		No peaks exceed the Max Peak Width threshold.
				One or more peaks exceed the Max Peak Width threshold.
CC	Control Concordance	Indicates if a positive, custom, or negative control produces the expected profile.  The expected profile is obtained from: <ul style="list-style-type: none"> <li>Positive Control – The Control Alleles in the panel used for analysis.</li> <li>Custom Control – The profile associated with the Profile ID in the Profile Manager.</li> <li>Negative Control – No peaks expected.</li> </ul> Applies only to Sample Type = Positive Control or Negative Control.  Weighting of the CC PQV is non-editable and set to 1.0.		Positive or Custom Control – Expected profile is detected.
				Negative Control – No peaks are detected.
				Positive or Custom Control – Expected profile is not detected.
				Negative Control – Peaks are detected.
			N/A	Sample Type = Sample or Allelic Ladder

Marker-level quality flags that are displayed in the Genotypes table after the analysis is complete (*continued*)




Acronym	Full Name	Function/Rule(s)	Flag Indicator	
SPK	Spike	<ul style="list-style-type: none"> <li>• <b>Allelic Ladders</b> – Indicates if spikes are detected within each extended marker range (no gaps are present between markers; the end point of each marker is extended past the marker definition in the panel to the beginning of the next marker).</li> <li>• <b>Samples</b> – Indicates if spikes are detected within a marker size range.</li> </ul> <p>The software uses a proprietary algorithm that detects spikes based on the peak morphology.</p>		No spikes are detected within the range.
				One or more spikes are detected within the range.
OVL	Overlapping Alleles	<p>Indicates if a labeled peak (allele or artifact) is within the size ranges of two neighboring markers.</p> <p>Applies only if you modify the default marker size ranges to overlap one another.</p>		No labeled peaks are detected within overlapping marker size ranges.
				One or more labeled peaks are detected within overlapping marker size ranges.

## Genotype Quality Assessment






For samples, the quality value system assigns the genotype quality (GQ) for each marker based on the individual marker quality flags.

For allelic ladders, the quality value system assigns a GQ for each marker based on the allelic ladder quality requirements, as described in “Allelic Ladder Quality Assessment” on page 45.

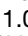









The GQ is used to determine the CGQ.

For samples, you can adjust the ranges that correspond to the Pass , Check , and Low Quality  ranges in the analysis method (SQ & GQ Settings).

Genotype Quality that is displayed in the Genotypes Table after the analysis is complete




Acronym	Full Name	Function/Rule(s)	Flag Indicator
GQ (samples)	Genotype Quality	<p>Indicates the genotype quality of the marker in the sample. The genotype quality for a sample marker is determined based on the presence of labeled peaks detected (after filtering) and the GQ weighting specified in the analysis method (SQ &amp; GQ Settings tab):</p> <ul style="list-style-type: none"> <li>If no labeled peaks are detected (and the sample is not a negative control), the GQ is set to 0.</li> <li>If one or more labeled peaks are detected, the GQ is initially set to 1 with a final value determined by the GQ weighting of individual marker-level quality values as specified in the analysis method using the formula:  <math display="block">GQ = 1 \times [(1 - \text{weight of flag 1}) \times (1 - \text{weight of flag 2}) \times (1 - \text{weight of flag } n)]</math>                     where:  <i>weight of flag 1</i> through <i>weight of flag n</i> represent the GQ weight for the following PQVs (applied only if the PQV is ): OS, BIN, PHR, LPH, AN, BD, CC, SPK, and OVL.</li> </ul>	 GQ is within the Passing range (Default = 0.75 to 1.0)
			 GQ is between the Passing range and Low Quality range (Default = 0.26 to 0.74)
			 GQ is within the Low Quality range (Default = 0.0 to 0.25) or the SQ for the sample is  .

Genotype Quality that is displayed in the Genotypes Table after the analysis is complete (*continued*)






Acronym	Full Name	Function/Rule(s)	Flag Indicator
GQ ( <i>continued</i> ) (samples)	Genotype Quality	<p><b>Weighting</b></p> <p>The weighting of individual flags affects the GQ. The CC weight is defined by the software at a value of 1.0. If CC is , GQ is set to 0 (  ). All other PQVs apply user-defined weights:</p> <ul style="list-style-type: none"> <li>• 1: If a PQV is triggered, it sets the GQ value to .</li> <li>• &gt; 0 &lt; 1: If a PQV is triggered, it reduces the GQ value according to the formula on the previous page. The higher the value, the greater the effect on GQ.</li> <li>• 0: No effect on the GQ calculation.</li> </ul> <p><b>Example</b></p> <p>A marker contains a  LPH and  OS flag, all other flags are . The GQ weight for LPH and OS are set to 0.3.</p> $GQ = 1 \times [(1 - 0.3) \times (1 - 0.3)]$ $GQ = 1 \times 0.49$ $GQ = 0.49$	
GQ (allelic ladders)	Genotype Quality	<p>Indicates the genotype quality of the marker in the allelic ladder. The genotype quality for an allelic ladder marker is determined using system-defined quality rules (described on <a href="#">page 45</a>) to ensure:</p> <ul style="list-style-type: none"> <li>• All expected peaks are present.</li> <li>• Peak height ratio of the first and second peak is greater than 50%.</li> <li>• No spikes are present in the extended marker range (within or between markers).</li> <li>• The peak height ratio between the lowest and highest peak is equal to or greater than 15%.</li> </ul> <p><b>IMPORTANT!</b> If the Allelic Ladder GQ Weighting for Spikes is set to 0 (off) (SQ &amp; GQ Settings tab in the analysis method), the GQ may be , even if spikes are present in the allelic ladder.</p>	<p> All allelic ladder quality requirements are met.</p> <p> One or more allelic ladder quality requirements are not met or the SQ for the sample is .</p>

## Sample-Level Quality Assessments












Sample-level quality assessment indicates the quality of the entire sample. Samples are grouped in the Analysis Summary based on the status of sample-level PQVs.

**Note:** Gray PQVs (  ) indicate that a marker in the sample was edited.

Sample-Level Quality Flags that are displayed in the Samples Table after the analysis is complete






Acronym	Full Name	Function/Rule(s)	Flag	
SOS	Sample Off-scale	Indicates if any fluorescence signal within the analysis range exceeds the detection threshold of the instrument.  <b>Note:</b> The detection limit varies on different instruments. The OS quality flag indicates if the CCD camera is saturated at any data point within the analysis range. It is not based on an absolute RFU value (peak height).		No fluorescence signal within the analysis range exceeds the detection limit of the instrument.
				One or more fluorescence signals within the analysis range exceed the detection limit of the instrument.
MIX	Mixed Source	Indicates a potential mixed source sample. A sample is considered a potential mixture if it meets either of the following conditions: <ul style="list-style-type: none"> <li>• <b>Condition 1:</b> Two or more markers contain three or more called alleles <i>or</i></li> <li>• <b>Condition 2:</b> <ul style="list-style-type: none"> <li>– One or more markers contain three or more called alleles <i>and</i></li> <li>– Another two or more markers have  PHR (peak height ratio, described on <a href="#">page 48</a>)</li> </ul> </li> </ul>		Neither condition is met.
				Either or both conditions are met, potential mixture.
			N/A	Sample Type = Allelic Ladder

Sample-Level Quality Flags that are displayed in the Samples Table after the analysis is complete (*continued*)

Acronym	Full Name	Function/Rule(s)	Flag	
OMR	Outside Marker Range	<p>Indicates if labeled peaks are detected between two marker size ranges defined in the panel. Does not indicate if labeled peaks are detected before the first marker or after the last marker.</p> <p><b>Note:</b> The default marker size ranges for neighboring markers do not overlap, but microvariants can fall between two marker size ranges.</p>		No labeled peaks are detected between marker size ranges.
				One or more labeled peaks are detected between marker size ranges.
SSPK	Sample Spike	<ul style="list-style-type: none"> <li>• <b>Allelic Ladders</b> – Indicates if spikes are detected within the sizing range.</li> <li>• <b>Samples</b> – Indicates if spikes are detected within or between two defined marker size ranges. Does not indicate if spikes are detected before the first marker or after the last marker.</li> </ul> <p>The software uses a proprietary algorithm that detects spikes based on the peak morphology.</p>		No spikes are detected within the range.
				One or more spikes are detected within the range.
CGQ (samples)	Composite Genotype Quality	Overall sample genotype quality indicator. Considers the individual marker GQ (Genotype Quality) values.		All individual marker GQ values are  .
				One or more individual marker GQ values are  .
				One or more individual marker GQ values are  .
				Manually overridden (see the GeneMapper® ID-X Software Help for information).



Sample-Level Quality Flags that are displayed in the Samples Table after the analysis is complete (*continued*)

Acronym	Full Name	Function/Rule(s)	Flag	
CGQ (allelic ladders)	Composite Genotype Quality	<p>Allelic ladder quality indicator. Considers the allelic ladder quality assessment (see <a href="#">page 45</a>). Other than SPK, does not consider individual marker GQ (Genotype Quality) values.</p> <p><b>Note:</b> Allelic ladder samples with  CGQ are not used to create bin offsets. For more information, see <a href="#">page 27</a>.</p> <p><b>IMPORTANT!</b> If the Allelic Ladder GQ Weighting for Spikes is set to 0 (off), or spike detection for the allelic ladder is disabled (SQ &amp; GQ Settings tab in the analysis method), the CGQ may be , even if spikes are present in the allelic ladder.</p>		All allelic ladder quality requirements are met for every marker.
				One or more allelic ladder quality requirements are not met for one or more markers.
				Manually overridden (see the GeneMapper® ID-X Software Help for information).





# Troubleshooting

A

Complete the following form before contacting Applied Biosystems Technical Support.

## Software Issues

Does the computer and operating system meet the required computer specifications?  
List any potential discrepancies.

Problem description:

Did you receive any error messages or alerts? If so, what did they say (if possible, take screen captures)? At what point did these errors/alerts occur?

Is the problem reproducible?



Steps taken prior to encountering or to reproduce the problem:

- 1.
- 2.
- 3.
- 4.
- 5.

---

# Bibliography

Baron, H., Fung, S., Aydin, A., Bahring, S., Luft, F.C., Schuster, H. 1996. Oligonucleotide ligation assay (OLA) for the diagnosis of familial hypercholesterolemia. *Nat Biotechnol.* 14(10):1279–82.

Brinkmann, B., Moller, A., and Wiegand, P. 1995. Structure of new mutations in 2 STR systems. *Int. J. Legal Med.* 107:201–203.

Gill, P., d'Aloja, E., Andersen, J., Dupuy, B., Jangblad, M., Johnsson, V., Kloosterman, A.D., Kratzer, A., Lareu, M.V., Meldegaard, M., Phillips, C., Pfitzinger, H., Rand, S., Sabatier, M., Scheithauer, R., Schmitter, H., Schneider, P., and Vide, M.C. 1997. Report of the European DNA Profiling Group (EDNAP): an investigation of the complex STR loci D21S11 and HUMFIBRA (FGA).

Grossman, P.D., Bloch, W., Brinson, E., Chang, C.C., Eggerding, F.A., Fung, S., Iovannisci, D.M., Woo, S., Win-Deen, E.S. 1994. High-density multiplex detection of nucleic acid sequences: oligonucleotide ligation assay and sequence-coded separation. *Nucleic Acids Res.* 22(21):4527–34.

Southern, E.M. 1979. Measurement of DNA length by gel electrophoresis. *Anal Biochem.* 100(2):319–23.



# Index

## Numerics

- 310, allelic ladder requirements 2
- 3100 and 3130 series, allelic ladder requirements 2

## A

- allele calling 34
- allele matching 26
- Allele Number 49
- alleles
  - naming convention 35
  - physical 28
  - that fall within virtual bins 28
- allelic ladder
  - analysis method for 3
  - averaging 29
  - CGQ 55
  - number per run, suggested 2
  - quality assessment 27, 45
  - reference bin sizes 27
  - requirements for accurate genotyping 2
  - sample type 3
- Amelogenin Cutoff 34
- AmpFISTR® Allelic Ladder and Binsets 27
- AN (allele number) 49
- analysis method for allelic ladders 3
- Analysis Requirement Not Met 42
- Analysis Requirements Summary 42
- analysis requirements, description 41
- ARNM (analysis requirements not met) 42
- assumptions for using this guide vii

## B

- baseline window 6

- BD (broad peak) 49
- BIN (out of bin allele) 47
- bin definitions 3
- bin offsetting 27, 29
- bold text, when to use vii
- Broad Peak 49

## C

- CC (control concordance) 49
- CGQ (composite genotype quality)
  - allelic ladder 55
  - samples and controls 54
- Control Concordance 49
- conventions
  - bold text vii
  - for describing menu commands viii
  - IMPORTANT! viii
  - in this guide vii
  - italic text vii
  - Notes viii
  - user attention words viii
- Cubic Spline Interpolation Method,
  - size-calling 24

## D

- disclaimer, license ii

## E

- examining peak definitions 5
- expert system, optimizing and validating 38

## F

- Find Missing Sample Files 41
- fluorescence signal threshold flag 53

## G

genotype PQVs, see marker PQVs

genotyping

allele calling 34

bin offsetting 27

overview 34

Global Cut-off value 34

Global Southern Method size-calling 21

Global Stutter 34

GQ (genotype quality)

description 51

displaying as number 39

sample 51, 52

weighting 52

## I

italic text, when to use vii

## L

Least Squares Method size-calling 22

license disclaimer ii

Local Southern Method size-calling 19

Low Peak Height 48

LPH (low peak height) 48

## M

marker margins, displaying offset 33

marker PQVs

AN (allele number) 49

BD (broad peak) 49

BIN (out of bin) 47

CC (control concordance) 49

GQ (genotype quality) 51, 52

LPH (low peak height) 48

MPH (max peak height) 48

OS (off-scale) 47

OVL (overlapping alleles) 50

PHR (peak height ratio) 48

SPK (spike) 50

marker-level quality assessments 47

marker-specific stutter ratios 34

Matrix Not Found 41

Max Peak Height 48

menu commands, conventions for describing viii

Minus A 34

MIX (mixture) 53

Mixed Source 53

MNF (matrix not found) 41

MPH (max peak height) 48

MSDSs, obtaining ix

## O

off-ladder alleles 35

Off-scale 47

offset bins 29

offset marker margins 33

OL allele calls 35

OMR (outside maker range) 54

optimizing peak detection sensitivity 11, 13

OS (off-scale) 47

Out of Bin 47

Outside Marker Range 54

Overlapping Alleles 50

OVL (overlap) 50

## P

peak definitions, examining 5

peak detection

baseline window 6

effects of extreme settings 13

optimizing sensitivity 11, 13

overview 4

parameters 4

peak window size 9

polynomial degree 9

sensitivity, optimizing 11, 13

slope threshold 13

slope thresholds peak start/end 13

smoothing 8

Peak Height Ratio 48

peak start, apex, and end marks 5

peak window size 9



PHR (peak height ratio) 48  
 polynomial degree  
   peak detection 9  
   varying 9  
   window size value 10  
 possible local sizing inaccuracy 25

## Q

quality assessments  
   allelic ladder 27  
   composite genotyping (CGQ) 36  
   genotype (GQ) 36  
   sizing (SQ) 25  
 quality value system  
   allelic ladder quality assessment 45  
   analysis requirements 41  
   checks and assessments 40  
   marker-level quality assessments 47  
   optimizing 38  
   overview 38  
   sample sizing quality 43  
   sample-level quality assessments 53

## R

Range Filter 34  
 ratio-matching 16  
 resolution, flag for loss of 49

## S

Sample File Not Found 41  
 Sample Off-scale 53  
 sample PQVs  
   CGQ (composite genotype quality) 54, 55  
   MIX (mixture) 53  
   OMR (outside marker range) 54  
   SOS (sample off-scale) 53  
   SQ (sizing quality) 43, 44  
   SSPK (sample spike) 54  
 Sample Spike 54  
 saturation flag 53  
 sensitivity, optimizing 11, 13

SFNF (sample file not found) 41  
 Show Peak Positions 5  
 size standard definition 16  
 Size Standard Not Found 41  
 size-calling  
   Cubic Spline Interpolation Method 24  
   Global Southern Method 21  
   Least Squares Method 22  
   Local Southern Method 19  
   overview 4  
 sizing  
   overview 16  
   quality (SQ). *See* SQ  
   size-calling curve and size-calling 18  
   size-matching 16  
   size-matching/size-calling algorithm 19, 21  
 Sizing Quality, *see* SQ  
 slope threshold  
   asymmetrical peak 15  
   peak end parameters 13  
   peak start parameters 13  
 smoothing 8  
 SOS (sample off-scale) 53  
 Spike 50  
 SPK (spike) 50  
 SQ (sizing quality)  
   description 43  
   displaying as number 39  
   weighting 43  
 SSNF (size standard not found) 41  
 SSPK (sample spike) 54  
 STRbase 28

## T

text conventions vii  
 training, information on ix  
 troubleshooting 57

## U

user attention words, described viii

## V

varying polynomial degree [9](#)

virtual bins [28](#)

## W

weighting

    GQ [52](#)

    SQ [43](#)

window size value, increasing [10](#)

### **Worldwide Sales and Support**

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches 150 countries on six continents. For sales office locations and technical support, please call our local office or refer to our Web site at [www.appliedbiosystems.com](http://www.appliedbiosystems.com).

Applied Biosystems is committed to providing the world's leading technology and information for life scientists.

### **Headquarters**

850 Lincoln Centre Drive  
Foster City, CA 94404 USA  
Phone: +1 650.638.5800  
Toll Free (In North America): +1 800.345.5224  
Fax: +1 650.638.5884

10/2009