

# AmpF $\ell$ STR<sup>™</sup> SGM Plus<sup>™</sup> PCR Amplification Kit

## USER GUIDE

Catalog Number 4307133

Publication Number 4309589

Revision J



**Manufacturer:** Thermo Fisher Scientific | 7 Kingsland Grange | Warrington, Cheshire WA1 4SR | United Kingdom

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Revision history:** Pub. No. 4309589

Revision	Date	Description
J	24 August 2018	Updated branding and trademarks, no technical changes.
H	August 2012	<ul style="list-style-type: none"><li>• Add validation experiments and results for buffer and enzyme kit component changes.</li><li>• Update instrument and software instructions.</li></ul>
G	March 2012	Change to copyright page information.
F	2011	Change to copyright page information.
E	2006	Change to copyright page information.
D	2006	—
C	2005	—
B	2001	—
A	1999	New document.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

**Trademarks:** All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license. AmpliTaq Gold is a trademark of Roche Molecular Systems, Inc. Windows and Windows Vista are trademarks of Microsoft Corporation. FTA is a trademark of Whatman International Limited. Minitab is a trademark of Minitab, Inc. Mac OS is a trademark of Apple, Inc.

©2018 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

<b>About This Guide</b> .....	<b>9</b>
Revision history .....	9
Purpose .....	9
User attention words .....	9
<b>■ CHAPTER 1 Overview</b> .....	<b>11</b>
Product overview .....	11
Purpose .....	11
Product description .....	11
About the primers .....	11
Loci amplified by the kit .....	11
Allelic ladder profile .....	12
Control DNA 007 profile .....	14
Workflow overview .....	15
Instrument and software overview .....	16
Data Collection and GeneMapper™ ID or ID-X Software .....	16
Instrument and software compatibility .....	16
About multicomponent analysis .....	16
How multicomponent analysis works .....	17
Materials and equipment .....	18
Kit contents and storage .....	18
Standards for samples .....	18
<b>■ CHAPTER 2 Perform PCR</b> .....	<b>19</b>
Required user-supplied reagents .....	19
DNA quantification .....	19
Importance of quantification .....	19
Methods of quantifying DNA .....	20
Prepare the amplification kit reactions .....	21
Select the correct PCR cycle number .....	22
Perform PCR .....	22
Amplification using bloodstained FTA™ cards .....	23

■	<b>CHAPTER 3 Electrophoresis</b>	<b>25</b>
	Allelic ladder requirements	25
	<b>Section 3.1 3100/3100-<i>Avant</i> and 3130/3130<i>xl</i> instruments</b>	<b>27</b>
	Set up the 3100/3100- <i>Avant</i> or 3130/3130 <i>xl</i> instrument for electrophoresis	27
	Reagents and parts	27
	Electrophoresis software setup and reference documents	27
	Prepare samples for electrophoresis on the 3100/3100- <i>Avant</i> or 3130/3130 <i>xl</i> instrument	28
	<b>Section 3.2 3500/3500xL Series instruments</b>	<b>29</b>
	Set up the 3500/3500xL instrument for electrophoresis	29
	Reagents and parts	29
	Electrophoresis software setup and reference documents	29
	Prepare samples for electrophoresis on the 3500/3500xL instrument	29
	<b>Section 3.3 310 Instrument</b>	<b>31</b>
	Set up the 310 instrument for electrophoresis	31
	Reagents and parts	31
	Electrophoresis software setup and reference documents	31
	Prepare samples for electrophoresis on the 310 instrument	31
■	<b>CHAPTER 4 Data Analysis</b>	<b>33</b>
	<b>Section 4.1 GeneMapper™ <i>ID</i> Software</b>	<b>33</b>
	Overview of GeneMapper™ <i>ID</i> Software	33
	Instruments	33
	Before you start	34
	Set up GeneMapper™ <i>ID</i> Software for data analysis	34
	File names	34
	Workflow	34
	Import panels and bins	35
	Create an analysis method	38
	General tab settings	39
	Allele tab settings	40
	Peak Detector tab settings	41
	Peak Quality tab settings	42
	Quality Flags tab settings	43
	Create size standard	43
	Analyze and edit sample files with GeneMapper™ <i>ID</i> Software	45
	Examine and edit a project	46
	For more information	46

<b>Section 4.2</b>	<b>GeneMapper™ <i>ID-X</i> Software</b>	<b>48</b>
Overview of GeneMapper™ <i>ID-X</i> Software		48
Instruments		48
Before you start		48
Set up GeneMapper™ <i>ID-X</i> Software for data analysis		49
Panel, bin, and stutter file version		49
Before using the software for the first time		49
Check panel, bin, and stutter file version		49
Import panels, bins, and marker stutter		50
Create an analysis method		55
General tab settings		56
Allele tab settings		57
Peak Detector tab settings		58
Peak Quality tab settings		59
SQ & GQ tab settings		60
Create size standard (optional)		60
Analyze and edit sample files with GeneMapper™ <i>ID-X</i> Software		62
Examine and edit a project		63
For more information		64
<b>■ CHAPTER 5</b>	<b>Experiments and Results</b>	<b>65</b>
<b>Section 5.1</b>	<b>Developmental Validation</b>	<b>66</b>
Overview		66
Importance of validation		66
Experiment conditions		66
Developmental validation		66
DAB 8.1.1 Developmental Validation		66
PCR components		66
Thermal cycler parameters		67
AmpliTaq Gold™ DNA Polymerase activation		67
Accuracy, reproducibility, and precision		67
DAB 8.1.2 Accuracy		67
Accuracy		68
Reproducibility		68
Precision and size windows		69
Extra peaks in the electropherogram		74
Overview		74
Stutter products		74
Addition of 3' A nucleotide		77
DAB 8.1.2.2 Mixture Studies		77

Characterization of loci .....	82
DAB 8.1.2.1 Documentation .....	82
Nature of the polymorphisms .....	82
Inheritance .....	83
Mapping .....	83
Population studies .....	83
Species specificity .....	83
DAB 8.1.2.2 Species Specificity .....	83
Nonhuman studies .....	84
Sensitivity .....	85
DAB 8.1.2.2 Sensitivity .....	85
Effect of DNA quantity on results .....	85
Stability .....	86
DAB 8.1.2.2 Stability .....	86
Lack of amplification of some loci .....	86
Differential and preferential amplification .....	87
Matrix studies .....	92
Population data .....	92
DAB 8.1.2.3 Population Data .....	92
DAB 8.1.2.3.1 Population Distribution Data .....	92
Overview .....	92
Population samples used in these studies .....	93
Allele frequencies .....	93
Mutation rate .....	99
Probability of identity .....	100
Probability of paternity exclusion .....	100
<b>Section 5.2 Performance Validation After Buffer and Enzyme Component</b>	
Replacement .....	102
Overview .....	102
Experiments .....	102
Reproducibility study .....	103
Intracolor balance .....	103
Stutter percentages .....	103
Artifacts .....	104
Sensitivity study .....	106
Mean peak height .....	106
DNA concentration and peak height .....	107
Allelic dropout .....	108
Genotype concordance .....	108
Inhibition study .....	108
Mean peak height, minimum mean peak height, and intracolor balance .....	108
Allelic dropout .....	112
Conclusions .....	112

■ <b>APPENDIX A</b>	<b>Troubleshooting</b>	<b>113</b>
■ <b>APPENDIX B</b>	<b>PCR Work Areas</b>	<b>115</b>
	Work area setup and lab design	115
	PCR setup work area	115
	Amplified DNA work area	116
■ <b>APPENDIX C</b>	<b>Ordering Information</b>	<b>117</b>
	Equipment and materials not included	117
■ <b>APPENDIX D</b>	<b>Safety</b>	<b>121</b>
	Chemical safety	122
	Biological hazard safety	123
	<b>Bibliography</b>	<b>125</b>
	<b>Documentation and Support</b>	<b>131</b>
	Related documentation	131
	Obtain SDSs	132
	Obtain support	132
	Limited Product Warranty	132
	<b>Index</b>	<b>133</b>



# About This Guide

---

**IMPORTANT!** Before using this product, read and understand the information the “Safety” appendix in this document.

---

## Revision history

Revision	Date	Description
A	1999	New document.
B	2001	—
C	2005	—
D	2006	—
E	2006	Change to limited licensing information.
F	2011	Change to limited licensing information.
G	March 2012	Change to copyright page information.
H	August 2012	<ul style="list-style-type: none"><li>• Add validation experiments and results for buffer and enzyme kit component changes.</li><li>• Update instrument and software instructions.</li></ul>

## Purpose

The Applied Biosystems *AmpFtSTR™ SGM Plus™ PCR Amplification Kit User Guide* provides information about the Applied Biosystems instruments, chemistries, and software associated with the AmpFtSTR™ SGM Plus™ PCR Amplification Kit.

## User attention words

Five user attention words may appear in this document. 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 or accurate chemistry kit use.

---



**CAUTION!** Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.

---



**WARNING!** Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.

---



**DANGER!** Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury.

---

- Product overview ..... 11
- Workflow overview ..... 15
- Instrument and software overview ..... 16
- Materials and equipment ..... 18

## Product overview

<b>Purpose</b>	The AmpF $\ell$ STR <sup>™</sup> SGM Plus <sup>™</sup> PCR Amplification Kit is a short tandem repeat (STR) multiplex assay that amplifies 10 tetranucleotide repeat loci (D3S1358, vWA, D16S539, D2S1338, D8S1179, D21S11, D18S51, D19S433, TH01, FGA) and the Amelogenin gender-determining marker in a single PCR amplification.
<b>Product description</b>	<p>The SGM Plus<sup>™</sup> Kit contains all the necessary reagents for the amplification of human genomic DNA.</p> <p>The reagents are designed for use with the following Applied Biosystems instruments:</p> <ul style="list-style-type: none"> <li>• Applied Biosystems 3100/3100-<i>Avant</i> Genetic Analyzer</li> <li>• Applied Biosystems 3130/3130<i>xl</i> Genetic Analyzer</li> <li>• Applied Biosystems 3500/3500<i>xL</i> Genetic Analyzer</li> <li>• Applied Biosystems 310 Genetic Analyzer</li> <li>• GeneAmp<sup>™</sup> PCR System 9700 with the Silver 96-Well Block</li> <li>• GeneAmp<sup>™</sup> PCR System 9700 with the Gold-plated Silver 96-Well Block</li> <li>• Veriti<sup>™</sup> 96-Well Thermal Cycler</li> </ul>
<b>About the primers</b>	The SGM Plus <sup>™</sup> Kit employs the same primer sequences for all loci common to other AmpF $\ell$ STR <sup>™</sup> kits (except the MiniFiler <sup>™</sup> kit), but does not include the degenerate unlabeled primer for the D8S1179 locus that was included in later AmpF $\ell$ STR <sup>™</sup> kits (for more information, see the <i>AmpF<math>\ell</math>STR<sup>™</sup> Identifiler PCR Amplification Kit User Guide</i> ).
<b>Loci amplified by the kit</b>	The following table shows the loci amplified, their chromosomal locations, and the corresponding fluorescent marker dyes. The AmpF $\ell$ STR <sup>™</sup> SGM Plus <sup>™</sup> Allelic Ladder is used to genotype the analyzed samples. The alleles contained in the allelic ladder and the genotype of the AmpF $\ell$ STR <sup>™</sup> Control DNA 007 are also listed in the table.

Table 1 SGM Plus™ Kit loci and alleles

Locus designation	Chromosome location	Alleles included in AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder	Dye label	Control DNA 007
D3S1358	3p	12, 13, 14, 15, 16, 17, 18, 19	5-FAM™	15, 16
vWA	12p12-pter	11, 12, 13, 14, 15, 16, 17, 18, 19, 20, 21, 22, 23, 24		14, 16
D16S539	16q24-qter	5, 8, 9, 10, 11, 12, 13, 14, 15		9, 10
D2S1338	2q35-37.1	15, 16, 17, 18, 19, 20, 21, 22, 23, 24, 25, 26, 27, 28		20, 23
Amelogenin	X: p22.1-22.3 Y: p11.2	X, Y	JOE™	X, Y
D8S1179†	8	8, 9, 10, 11, 12, 13, 14, 15, 16, 17, 18, 19		12, 13
D21S11	21q11.2-q21	24, 24.2, 25, 26, 27, 28, 28.2, 29, 29.2, 30, 30.2, 31, 31.2, 32, 32.2, 33, 33.2, 34, 34.2, 35, 35.2, 36, 37, 38		28, 31
D18S51	18q21.3	7, 9, 10, 10.2, 11, 12, 13, 13.2, 14, 14.2, 15, 16, 17, 18, 19, 20, 21, 22, 23, 24, 25, 26, 27		12, 15
D19S433	19q12-13.1	9, 10, 11, 12, 12.2, 13, 13.2, 14, 14.2, 15, 15.2, 16, 16.2, 17, 17.2	NED™	14, 15
TH01	11p15.5	4, 5, 6, 7, 8, 9, 9.3, 10, 11, 13.3		7, 9.3
FGA	4q28	17, 18, 19, 20, 21, 22, 23, 24, 25, 26, 26.2, 27, 28, 29, 30, 30.2, 31.2, 32.2, 33.2, 42.2, 43.2, 44.2, 45.2, 46.2, 47.2, 48.2, 50.2, 51.2		24, 26

† In some literature references, this locus is designated as D6S502.

## Allelic ladder profile

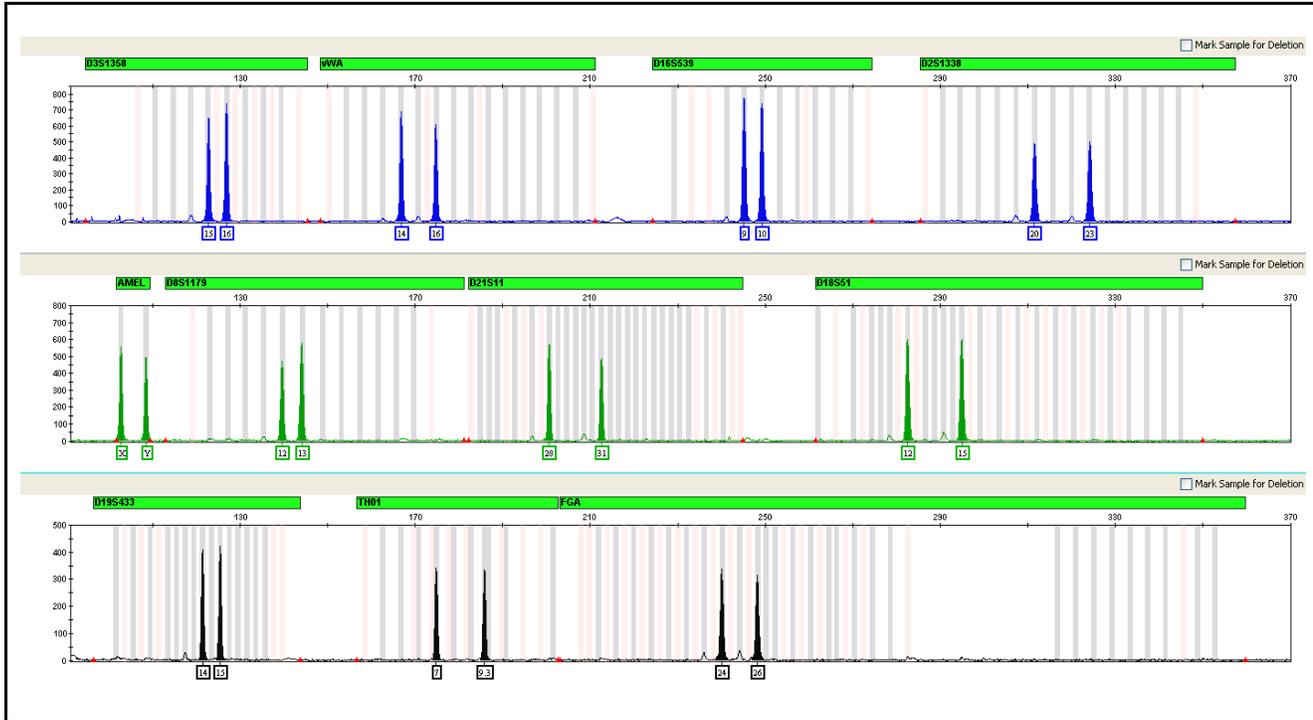
Figure 1 shows the allelic ladder for the SGM Plus™ Kit. See “Allelic ladder requirements” on page 25 for information on ensuring accurate genotyping.



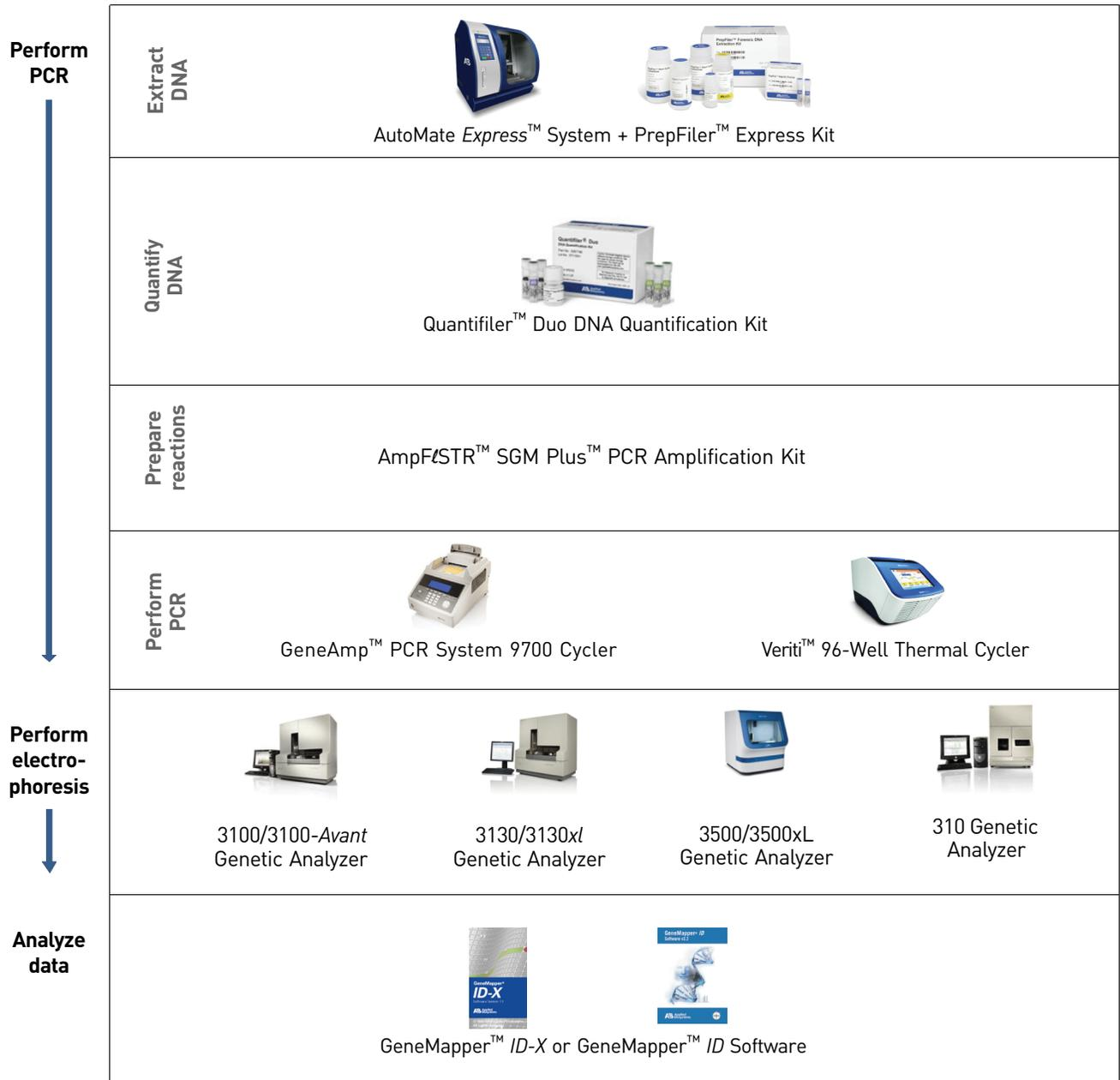
### Control DNA 007 profile

Figure 2 shows amplification of Control DNA 007 using the SGM Plus™ Kit.

Figure 2 1 ng of Control DNA 007 amplified with the SGM Plus™ Kit and analyzed on the Applied Biosystems 3130xl Genetic Analyzer



# Workflow overview



## Instrument and software overview

This section provides information about the Data Collection Software versions required to run the AmpF $\lambda$ STR™ SGM Plus™ PCR Amplification Kit on specific instruments.

### Data Collection and GeneMapper™ ID or ID-X Software

The Data Collection Software provides instructions to firmware running on the instrument and displays instrument status and raw data in real time. As the instrument measures sample fluorescence with its detection system, the Data Collection Software collects the data and stores it. The Data Collection Software stores information about each sample in a sample file (.fsa), which is then analyzed by the GeneMapper™ ID or ID-X Software.

### Instrument and software compatibility

**Table 2** Software specific to each instrument

Instrument	Operating system	Data Collection Software	Analysis software
3500/3500xL	<ul style="list-style-type: none"> <li>Windows™ XP</li> <li>Windows Vista™</li> </ul>	3500 Series Data Collection Software v1.0	GeneMapper™ ID-X Software v1.2 or higher
3130/3130xl	Windows™ XP	3.0	<ul style="list-style-type: none"> <li>GeneMapper™ ID Software v3.2.1</li> <li>and</li> <li>GeneMapper™ ID-X Software v1.0.1 or higher</li> </ul>
3100/3100-Avant	Windows™ NT	1.1 (3100) 1.0 (3100-Avant)	
	Windows™ 2000	2.0	
310	Windows™ XP	3.1	
	<ul style="list-style-type: none"> <li>Windows™ NT</li> <li>Windows™ 2000</li> </ul>	3.0	

**Note:** We conducted validation studies for the AmpF $\lambda$ STR™ SGM Plus™ PCR Amplification Kit using the Applied Biosystems 310 Genetic Analyzer running Mac Operating System™. This configuration is now obsolete.

### About multicomponent analysis

Applied Biosystems fluorescent multi-color dye technology allows the analysis of multiple loci, including loci that have alleles with overlapping size ranges. Alleles for overlapping loci are distinguished by labeling locus-specific primers with different colored dyes.

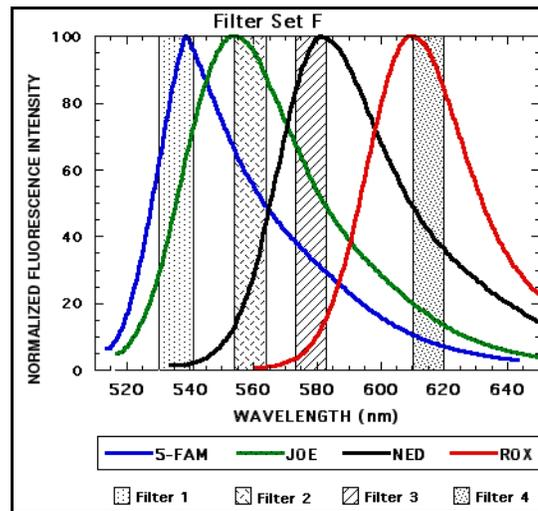
Multicomponent analysis is the process that separates the four different fluorescent dye colors into distinct spectral components. The three dyes used in the SGM Plus™ Kit to label samples are 5-FAM™, JOE™, and NED™ dyes. The fourth dye, ROX™ dye, is used to label the GeneScan™ 500 ROX™ Size Standard.

## How multicomponent analysis works

Each of these fluorescent dyes emits its maximum fluorescence at a different wavelength. During data collection on the Applied Biosystems and Applied Biosystems instruments, the fluorescence signals are separated by diffraction grating according to their wavelengths and projected onto a charge-coupled device (CCD) camera in a predictably spaced pattern. The 5-FAM™ dye emits at the shortest wavelength and it is displayed as blue, followed by JOE™ (green), NED™ (yellow), and ROX™ (red) dyes.

Although each of these dyes emits its maximum fluorescence at a different wavelength, there is some overlap in the emission spectra between the dyes (Figure 3). The goal of multicomponent analysis is to correct for spectral overlap.

Figure 3 Emission spectra of the four dyes used in the AmpFtSTR™ SGM Plus™ Kit



## Materials and equipment

### Kit contents and storage

The AmpF $\phi$ STR™ SGM Plus™ PCR Amplification Kit (Part no. 4307133) contains materials sufficient to perform 100 amplifications at 50  $\mu$ L/amplification.

**IMPORTANT!** The fluorescent dyes attached to the primers are light sensitive. Protect the primer set, amplified DNA, allelic ladder, and size standard from light when not in use. Keep freeze-thaw cycles to a minimum.

Component	Description	Volume	Storage
AmpF $\phi$ STR™ PCR Reaction Mix	Contains MgCl <sub>2</sub> , dNTPs, and bovine serum albumin in buffer with 0.05% sodium azide.	2 tubes, 1.1 mL each	-15 to -25°C on receipt, 2 to 8°C after initial use
AmpF $\phi$ STR™ SGM Plus™ Primer Set	Contains fluorescently labeled primers and non-labeled primers.	1 tube, 1.1 mL	
AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder	Contains amplified alleles. See <a href="#">Table 1 on page 12</a> for a list of alleles included in the allelic ladder.	1 tube, 0.05 mL	
AmpF $\phi$ STR™ Control DNA 007	Contains 0.10 ng/ $\mu$ L human male 007 DNA in 0.05% sodium azide and buffer†. See <a href="#">Table 1 on page 12</a> for profile.	1 tube, 0.3 mL	
AmpliQ Gold™ DNA Polymerase	Contains enzyme, with an activity of 5 U/ $\mu$ L.	2 tubes, 0.05 mL/tube	-15 to -25°C

† The AmpF $\phi$ STR™ Control DNA 007 is included at a concentration appropriate to its intended use as an amplification control (i.e., to provide confirmation of the capability of the kit reagents to generate a profile of expected genotype). The AmpF $\phi$ STR™ Control DNA 007 is not designed to be used as a DNA quantitation control, and you may see variation from the labelled concentration when quantitating aliquots of the AmpF $\phi$ STR™ Control DNA 007.

### Standards for samples

For the SGM Plus™ Kit, the panel of standards needed for PCR amplification, PCR product sizing, and genotyping are:

- **AmpF $\phi$ STR™ Control DNA 007** – A positive control for evaluating the efficiency of the amplification step and STR genotyping using the AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder.
- **GeneScan™ 500 ROX™ Size Standard** – Used for obtaining sizing results. This standard, which has been evaluated as an internal size standard, yields precise sizing results for SGM Plus™ Kit PCR products. Order the GeneScan™ 500 ROX™ Size Standard (Part no. 4322682) separately.
- **AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder** – Allelic ladder developed by Life Technologies for accurate characterization of the alleles amplified by the SGM Plus™ Kit. The AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder contains most of the alleles reported for the 10 autosomal loci. Refer to [Table 1 on page 12](#) for a list of the alleles included in the AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder.

■ Required user-supplied reagents . . . . .	19
■ DNA quantification . . . . .	19
■ Prepare the amplification kit reactions . . . . .	21
■ Select the correct PCR cycle number . . . . .	22
■ Perform PCR . . . . .	22
■ Amplification using bloodstained FTA™ cards . . . . .	23

## Required user-supplied reagents

In addition to the SGM Plus™ Kit reagents, the use of low-TE buffer (10 mM Tris, 0.1 mM EDTA, pH 8.0) is recommended. You can prepare the buffer as described in the procedure below or order it from Teknova (Cat # T0223).

To prepare low-TE buffer:

- Mix together:
  - 10 mL of 1 M Tris-HCl, pH 8.0
  - 0.2 mL of 0.5 M EDTA, pH 8.0
  - 990 mL glass-distilled or deionized water

**Note:** Adjust the volumes accordingly for specific needs.
- Aliquot and autoclave the solutions.
- Store at room temperature.

## DNA quantification

### Importance of quantification

Quantifying the amount of DNA in a sample before amplification allows you to determine whether or not sufficient DNA is present to permit amplification and to calculate the optimum amount of DNA to add to the reaction. The optimum amount of DNA for the SGM Plus™ Kit is 1.0–2.5 ng in a maximum input volume of 20 µL for 28 PCR cycles.

If too much DNA is added to the PCR reaction, then the increased amount of PCR product that is generated can result in:

- Fluorescence intensity that exceeds the linear dynamic range for detection by the instrument (“off-scale” data). Off-scale data are problematic because:
  - Quantitation (peak height and area) for off-scale peaks is not accurate. For example, an allele peak that is off-scale can cause the corresponding stutter peak to appear higher in relative intensity, thus increasing the calculated percent stutter.
  - Multicomponent analysis of off-scale data is not accurate, and it results in poor spectral separation (“pull-up”).
- Incomplete A-nucleotide addition.

When the total number of allele copies added to the PCR is extremely low, allelic dropout can occur resulting in a partial profile.

## Methods of quantifying DNA

Life Technologies provides several kits for quantifying DNA in samples. See the reference cited in the following table for details about these kits.

Product	Description
Quantifiler™ Human DNA Quantification Kit (Part no. 4343895)  <i>and</i> Quantifiler™ Y Human Male DNA Quantification Kit (Part no. 4343906)  For more information, see <i>Quantifiler™ Human DNA Quantification Kits User's Manual</i> (Pub. no. 4344790)	<p><b>Properties:</b></p> <p>The Quantifiler™ Human and Quantifiler™ Y Human Male Kits are highly specific for human DNA, and they individually detect total human or male DNA, respectively. The kits detect single-stranded and degraded DNA.</p> <p><b>How they work:</b></p> <p>The Quantifiler™ DNA Quantification Kits consist of target-specific and internal control 5' nuclease assays.</p> <p>The Quantifiler™ Human and Quantifiler™ Y Human Male Kits contain different target-specific assays (human DNA or human male DNA, respectively) that each consist of two locus-specific PCR primers and one TaqMan™ MGB probe labeled with FAM™ dye for detecting the amplified sequence. The kits each contain a separate internal PCR control (IPC) assay, which consists of an IPC template DNA (a synthetic sequence not found in nature), two primers for amplifying the IPC template, and one TaqMan™ MGB probe labeled with VIC™ dye for detecting the amplified IPC.</p>
Quantifiler™ Duo DNA Quantification Kit (Part no. 4387746)  For more information, see <i>Quantifiler™ Duo DNA Quantification Kit User's Manual</i> (Pub. no. 4391294)	<p><b>Properties:</b></p> <p>The Quantifiler™ Duo Kit is highly specific for human DNA. This kit combines the detection of both total human and male DNA in one PCR reaction. The kit detects single-stranded and degraded DNA.</p> <p><b>How it works:</b></p> <p>The Quantifiler™ Duo DNA Quantification Kit consists of target-specific and internal control 5' nuclease assays.</p> <p>The Quantifiler™ Duo kit combines two human-specific assays in one PCR reaction (for total human DNA and human male DNA). The two human DNA specific assays each consist of two PCR primers and a TaqMan™ probe. The TaqMan™ probes for the human DNA and human male DNA assays are labeled with VIC™ and FAM™ dyes, respectively. In addition, the kit contains an internal PCR control (IPC) assay similar in principle to that used in the other Quantifiler kits, but labeled with NED™ dye.</p>

## Prepare the amplification kit reactions

1. Calculate the volume of each component needed to prepare the reactions, using the table below.

DNA sample	Volume per reaction
AmpF $\lambda$ STR™ PCR Reaction Mix	21.0 $\mu$ L
AmpF $\lambda$ STR™ SGM Plus™ Primer Set	11.0 $\mu$ L
AmpliTaq Gold™ DNA Polymerase	1.0 $\mu$ L

**Note:** The volumes above include a slight overfill to provide excess volume for the loss that occurs during reagent transfers.

2. Prepare reagents. Thaw the AmpF $\lambda$ STR™ PCR Reaction Mix and the SGM Plus™ Primer Set, then vortex all reagent tubes including the enzyme for 3 seconds and centrifuge briefly before opening the tubes.

**IMPORTANT!** Thawing is required only during first use of the kit. After first use, reagents are stored at 2 to 8°C and, therefore, they do not require subsequent thawing. Do not refreeze these reagents.

3. Pipette the required volumes of components into an appropriately sized polypropylene tube to create a master mix.
4. Vortex the master mix for 3 seconds, then centrifuge briefly.
5. Dispense 30  $\mu$ L of the reaction mix into each reaction well of a MicroAmp™ Optical 96-Well Reaction Plate or each MicroAmp™ tube.
6. Prepare the DNA samples:

DNA sample	To prepare...
Negative control	Add 20 $\mu$ L of low-TE buffer (10mM Tris, 0.1mM EDTA, pH 8.0).
Test sample	Dilute a portion of the test DNA sample with low-TE buffer so that 1.0–2.5 ng of total DNA is in a final volume of 20 $\mu$ L. Add 20 $\mu$ L of the diluted sample to the reaction mix.
Positive control	Add 20 $\mu$ L of 007 control DNA (0.1 ng/ $\mu$ L).

The final reaction volume (sample or control plus master mix) is 50  $\mu$ L.

7. Seal the MicroAmp™ Optical 96-Well Reaction Plate with MicroAmp™ Clear Adhesive Film or MicroAmp™ Optical Adhesive Film, or cap the tubes.
8. Centrifuge the tubes or plate at 3000 rpm for ~20 seconds in a tabletop centrifuge (with plate holders if using 96-well plates).
9. Amplify the samples in a GeneAmp™ PCR System 9700 with the silver or gold-plated silver 96-well block or a Veriti™ 96-Well Thermal Cycler.

**Note:** The SGM Plus™ Kit is not validated for use with the GeneAmp PCR System 9700 with the aluminium 96-well block. Use of this thermal cycling platform may adversely affect performance of the SGM Plus™ Kit.

## Select the correct PCR cycle number

All AmpF $\Delta$ STR™ kits are optimized for a specific number of amplification cycles to deliver well-balanced and high quality results. However, increases in the number of low-level DNA samples being submitted for analysis have prompted many laboratories to evaluate increasing the number of amplification cycles to increase the sensitivity of the assay. Before increasing the cycle number, perform a comprehensive validation study to establish new performance criteria for the higher cycle number.

Higher cycle numbers can cause the following to occur:

- Exaggerated stochastic effects resulting from low DNA input amounts
- Greater difference between the presence and absence of an allele
- Greater heterozygote peak imbalance
- Possible differences in expected stutter position and percentage
- Possible increase in artifacts and/or background in the profile to accompany the increase in sample allele signal

The AmpF $\Delta$ STR™ SGM Plus™ PCR Amplification Kit is optimized for 28 cycles of amplification only.

The results of developmental validation studies are shown in [“Developmental Validation” on page 66](#).

## Perform PCR

1. Program the thermal cycling conditions:
  - When using the GeneAmp™ PCR System 9700 with either 96-well silver or gold-plated silver block, select the **9600 Emulation Mode**.
  - When using the Veriti™ 96-Well Thermal Cycler, refer to the following document for instructions on how to configure the Veriti instrument to run in the 9600 Emulation Mode: *User Bulletin: Veriti™ 96-Well Thermal Cycler AmpF $\Delta$ STR™ Kit Validation* (Pub. no.4440754).

Initial incubation step	Melt	Anneal	Extend	Final extension	Final hold
HOLD	CYCLE (28)			HOLD	HOLD
95°C 11 min	94°C 1 min	59°C 1 min	72°C 1 min	60°C 45 min	25°C ∞

2. Load the plate into the thermal cycler and close the heated cover.

---

**IMPORTANT!** If using the 9700 thermal cycler with silver or gold-plated silver block and adhesive clear film instead of caps to seal the plate wells, be sure to place a MicroAmp™ compression pad (Part no. 4312639) on top of the plate to prevent evaporation during thermal cycling. The Veriti™ Thermal Cycler does not require a compression pad.

---

3. Start the run.

- On completion of the run, store the amplified DNA and protect from light.

If you are storing the DNA...	Then place at...
< 2 weeks	2 to 8°C
> 2 weeks	-15 to -25°C

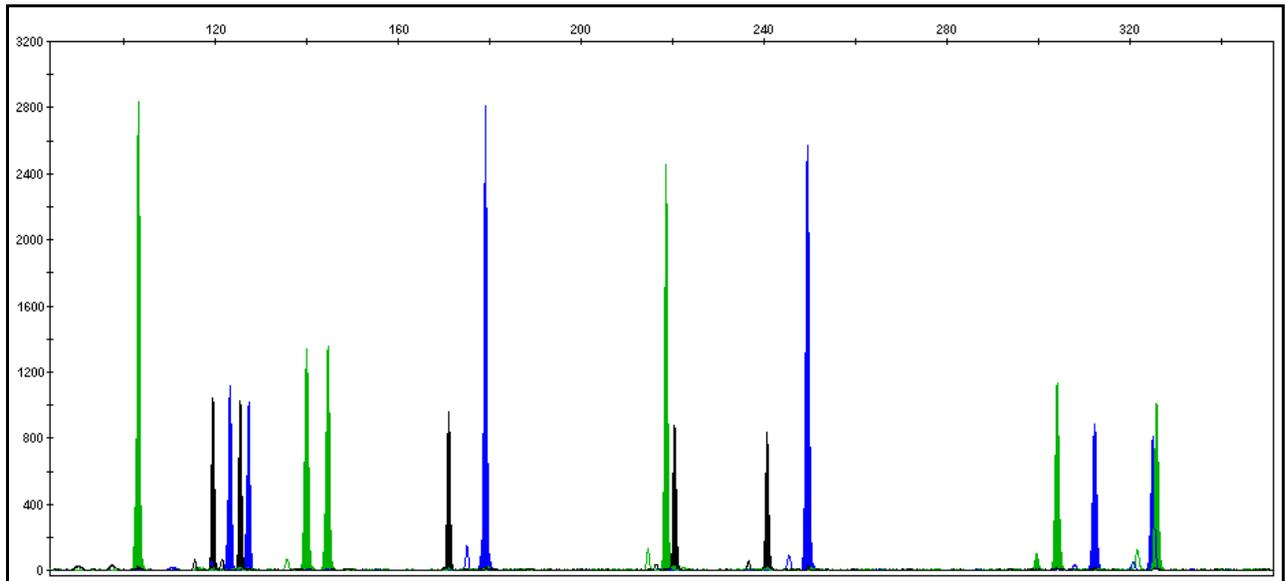
**IMPORTANT!** Store the amplified products so that they are protected from light.

## Amplification using bloodstained FTA™ cards

FTA™ cards can be useful for collecting, storing, and processing biological samples. A small punch disc of the card containing the sample can be placed directly into an amplification tube, purified, and amplified, without transferring the disc. Our studies indicate that a 1.2-mm bloodstained disc contains approximately 5–20 ng DNA. An appropriate cycle number for this high quantity of DNA is 25 cycles as determined by our validation studies. However, it is recommended that each laboratory determine the optimum cycle number based on internal validation studies.

In the example shown in [Figure 4](#), a 1.2-mm disc of a bloodstained FTA card was purified using three washes with FTA Purification Reagent and two washes with 1× low-TE buffer. The purified punch disc was then amplified in a MicroAmp™ tube for 25 cycles.

**Figure 4** AmpFtSTR™ SGM Plus™ Kit PCR Amplification Kit results from a 1.2-mm FTA bloodstain disc (25-cycle amplification), analyzed on the Applied Biosystems 3130xl Genetic Analyzer





■ Allelic ladder requirements . . . . .	25
■ 3100/3100-Avant and 3130/3130xl instruments . . . . .	27
Set up the 3100/3100-Avant or 3130/3130xl instrument for electrophoresis . . . .	27
Prepare samples for electrophoresis on the 3100/3100-Avant or 3130/3130xl instrument . . . . .	28
■ 3500/3500xL Series instruments . . . . .	29
Set up the 3500/3500xL instrument for electrophoresis . . . . .	29
Prepare samples for electrophoresis on the 3500/3500xL instrument. . . . .	29
■ 310 Instrument . . . . .	31
Set up the 310 instrument for electrophoresis . . . . .	31
Prepare samples for electrophoresis on the 310 instrument . . . . .	31

## Allelic ladder requirements

To accurately genotype samples, you must run an allelic ladder sample along with the unknown samples.

Instrument	Number of allelic ladders to run	One injection equals	Number of samples per allelic ladder(s)
3100-Avant or 3130	1 per 4 injections	4 samples	15 samples + 1 allelic ladder
3100 or 3130xl	1 per injection	16 samples	15 samples + 1 allelic ladder
3500	1 per 3 injections	8 samples	23 samples + 1 allelic ladder
3500xL	1 per injection	24 samples	23 samples + 1 allelic ladder
310	1 per 10 injections	1 sample	9 samples + 1 allelic ladder

**IMPORTANT!** Variation in laboratory temperature can cause changes in fragment migration speed and sizing variation between both single- and multiple-capillary runs (with larger size variations seen between samples injected in multiple-capillary runs). We recommend the above frequency of allelic ladder injections, which should account for normal variation in run 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 run under the same conditions as the samples because size values obtained for the same sample can differ between instrument platforms because of different polymer matrices and electrophoretic conditions.

## Section 3.1 3100/3100-Avant and 3130/3130xl instruments

### Set up the 3100/3100-Avant or 3130/3130xl instrument for electrophoresis

**Reagents and parts** “Ordering Information” on page 117 lists the required materials not supplied with the AmpFtSTR™ SGM Plus™ PCR Amplification Kit.

**IMPORTANT!** The fluorescent dyes attached to the primers are light sensitive. Protect the primer set, amplified DNA, allelic ladder, and size standard from light when not in use. Keep freeze-thaw cycles to a minimum.

#### Electrophoresis software setup and reference documents

The following table lists Data Collection Software and the run modules that can be used to analyze SGM Plus™ Kit PCR products. For details on the procedures, refer to the documents listed in the table.

Genetic Analyzer	Data Collection Software	Operating System	Run modules and conditions	References
Applied Biosystems 3130/3130xl	3.0	Windows™ XP	<ul style="list-style-type: none"> <li>HIDFragmentAnalysis36_POP4_1</li> <li>Injection conditions:                             <ul style="list-style-type: none"> <li>3130 = 3 kV/5 sec</li> <li>3130xl = 3 kV/10 sec</li> </ul> </li> <li>Dye Set F</li> </ul>	<i>Applied Biosystems 3130/3130xl Genetic Analyzers Using Data Collection Software v3.0, Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin</i> (Pub. no. 4363787)
Applied Biosystems 3100	2.0	Windows™ 2000	<ul style="list-style-type: none"> <li>HIDFragmentAnalysis36_POP4_1</li> <li>Injection condition: 3kV/10 sec</li> <li>Dye Set F</li> </ul>	<i>3100/3100-Avant Genetic Analyzers Using Data Collection Software v2.0, Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin</i> (Pub. no. 4350218)
	1.1	Windows™ NT	<ul style="list-style-type: none"> <li>GeneScan36_POP4DyeSetF</li> <li>Injection condition: 3kV/10 sec</li> <li>GS500Analysis.gsp</li> </ul>	<i>3100/3100-Avant Genetic Analyzers Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin</i> (Pub. no. 4332345)
Applied Biosystems 3100-Avant	1.0	Windows™ NT	<ul style="list-style-type: none"> <li>GeneScan36A_POP4DyeSetFModule</li> <li>Injection condition: 3 kV/5sec</li> <li>GS500Analysis.gsp</li> </ul>	<i>3100/3100-Avant Genetic Analyzers Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin</i> (Pub. no. 4332345)

## Prepare samples for electrophoresis on the 3100/3100-Avant or 3130/3130xl instrument

Prepare the samples for electrophoresis on the 3100/3100-Avant or 3130/3130xl instrument immediately before loading.

1. Calculate the volume of Hi-Di™ Formamide and size standard needed to prepare the samples:

Reagent	Volume per reaction
GeneScan™ 500 ROX™ Size Standard	0.5 µL
Hi-Di™ Formamide	8.5 µL

**Note:** Include additional samples in your calculations to provide excess volume for the loss that occurs during reagent transfers.

---

**IMPORTANT!** The volume of size standard indicated in the table is a suggested amount. Determine the appropriate amount of size standard based on your experiments and results.

---

2. Pipette the required volumes of components into an appropriately sized polypropylene tube.
3. Vortex the tube, then centrifuge briefly.
4. Into each well of a MicroAmp™ Optical 96-Well Reaction Plate, add:
  - 9 µL of the formamide:size standard mixture
  - 1 µL of PCR product or allelic ladder

**Note:** For blank wells, add 10 µL of Hi-Di™ Formamide.
5. Seal the reaction plate with appropriate septa, then centrifuge the plate to ensure that the contents of each well are collected at the bottom.
6. Heat the reaction plate in a thermal cycler for 3 minutes at 95°C.
7. Immediately place the plate on ice for 3 minutes.
8. Prepare the plate assembly, then place on the autosampler.
9. Ensure that a plate record is completed and link the plate record to the plate.
10. Start the electrophoresis run.

## Section 3.2 3500/3500xL Series instruments

### Set up the 3500/3500xL instrument for electrophoresis

**Reagents and parts** “Ordering Information” on page 117 lists the required materials not supplied with the AmpF<sub>STR</sub><sup>™</sup> SGM Plus<sup>™</sup> PCR Amplification Kit.

**IMPORTANT!** The fluorescent dyes attached to the primers are light sensitive. Protect the primer set, amplified DNA, allelic ladder, and size standard from light when not in use. Keep freeze-thaw cycles to a minimum.

#### Electrophoresis software setup and reference documents

The following table lists Data Collection Software and the run modules that can be used to analyze SGM Plus<sup>™</sup> Kit PCR products. For details on the procedures, refer to the documents listed in the table.

Genetic Analyzer	Data Collection Software	Operating System	Run modules and conditions	References
Applied Biosystems 3500	3500 Data Collection Software v1.0	Windows <sup>™</sup> XP	<ul style="list-style-type: none"> <li>HID36_POP4</li> <li>Injection conditions: 1.2kV/15 sec</li> <li>Dye Set F</li> </ul>	<i>Applied Biosystems 3500/3500xL Genetic Analyzer User Guide</i> (Pub. no. 4401661)
Applied Biosystems 3500xL		Windows Vista <sup>™</sup>	<ul style="list-style-type: none"> <li>HID36_POP4</li> <li>Injection conditions: 1.2kV/24 sec</li> <li>Dye Set F</li> </ul>	<i>Applied Biosystems 3500 and 3500xL Genetic Analyzers Quick Reference Card</i> (Pub. no. 4401662)

### Prepare samples for electrophoresis on the 3500/3500xL instrument

Prepare the samples for electrophoresis on the 3500/3500xL instrument immediately before loading.

1. Calculate the volume of Hi-Di<sup>™</sup> Formamide and size standard needed to prepare the samples:

Reagent	Volume per reaction
GeneScan <sup>™</sup> 500 ROX <sup>™</sup> Size Standard	0.5 µL
Hi-Di <sup>™</sup> Formamide	8.5 µL

**Note:** Include additional samples in your calculations to provide excess volume for the loss that occurs during reagent transfers.

---

**IMPORTANT!** The volume of size standard indicated in the table is a suggested amount. Determine the appropriate amount of size standard based on your results and experiments.

---

2. Pipette the required volumes of components into an appropriately sized polypropylene tube.
3. Vortex the tube, then centrifuge briefly.
4. Into each well of a MicroAmp™ Optical 96-Well Reaction Plate, or each MicroAmp™ optical strip tube, add:
  - 9 µL of the formamide:size standard mixture
  - 1 µL of PCR product or allelic ladder

**Note:** For blank wells, add 10 µL of Hi-Di™ Formamide.
5. Seal the reaction plate or strip tubes with the appropriate septa, then centrifuge to ensure that the contents of each well are collected at the bottom.
6. Heat the reaction plate or strip tubes in a thermal cycler for 3 minutes at 95°C.
7. Immediately put the plate or strip tubes on ice for 3 minutes.
8. Prepare the plate assembly, then place on the autosampler.
9. Ensure that a plate record is completed and link the plate record to the plate.
10. Start the electrophoresis run.

## Section 3.3 310 Instrument

### Set up the 310 instrument for electrophoresis

**Reagents and parts** “Ordering Information” on page 117 lists the required materials not supplied with the AmpF $\ell$ STR™ SGM Plus™ PCR Amplification Kit.

---

**IMPORTANT!** The fluorescent dyes attached to the primers are light sensitive. Protect the primer set, amplified DNA, allelic ladder, and size standard from light when not in use. Keep freeze-thaw cycles to a minimum.

---

#### Electrophoresis software setup and reference documents

The following table lists Data Collection Software and the run modules that can be used to analyze SGM Plus™ Kit PCR products. For details on the procedures, refer to the documents listed in the table.

Data Collection Software	Operating System	Run modules and conditions	References
3.1 <sup>†</sup> or 3.0 <sup>†</sup>	Windows™ XP or Windows™ NT and Windows™ 2000	<ul style="list-style-type: none"> <li>GS STR POP4 (1mL)</li> </ul> Injection condition: 15 kV/5 sec	<i>310 Genetic Analyzer User's Manual (Windows)</i> [Pub. no. 4317588]  <i>310 Protocols for Processing AmpF<math>\ell</math>STR™ PCR Amplification Kit Products with Microsoft Windows NT Operating System: User Bulletin</i> [Pub. no. 4341742]

<sup>†</sup> We conducted concordance studies for the SGM Plus™ Kit using this configuration.

### Prepare samples for electrophoresis on the 310 instrument

Prepare the samples for electrophoresis on the 310 instrument immediately before loading.

1. Calculate the volume of Hi-Di™ Formamide and size standard needed to prepare the samples:

Reagent	Volume per reaction
GeneScan™ 500 ROX™ Size Standard	0.75 $\mu$ L
Hi-Di™ Formamide	24.5 $\mu$ L

**Note:** Include additional samples in your calculations to provide excess volume for the loss that occurs during reagent transfers.

---

**IMPORTANT!** The volume of size standard indicated in the table is a suggested amount. Determine the appropriate amount of size standard based on your results and experiments.

---

2. Pipette the required volumes of components into an appropriately sized polypropylene tube.
3. Vortex the tube, then centrifuge briefly.
4. Into each 0.2 mL or 0.5 mL sample tube, add:
  - 25  $\mu$ L of the formamide: size standard mixture
  - 1.5  $\mu$ L of PCR product or allelic ladder

**Note:** For blank wells, add 25  $\mu$ L of Hi-Di™ Formamide.
5. Seal the tubes with the appropriate septa, then briefly centrifuge to ensure that the contents of each tube are mixed and collected at the bottom.
6. Heat the tubes in a thermal cycler for 3 minutes at 95°C.
7. Immediately place the tubes on ice for 3 minutes.
8. Place the sample tray on the autosampler.
9. Ensure that an injection list is prepared.
10. Start the electrophoresis run.

■ GeneMapper™ ID Software .....	33
Overview of GeneMapper™ ID Software .....	33
Set up GeneMapper™ ID Software for data analysis .....	34
Analyze and edit sample files with GeneMapper™ ID Software .....	45
Examine and edit a project .....	46
For more information .....	46
■ GeneMapper™ ID-X Software .....	48
Overview of GeneMapper™ ID-X Software .....	48
Set up GeneMapper™ ID-X Software for data analysis .....	49
Analyze and edit sample files with GeneMapper™ ID-X Software .....	62
Examine and edit a project .....	63
Examine and edit a project .....	63

## Section 4.1 GeneMapper™ ID Software

### Overview of GeneMapper™ ID Software

GeneMapper™ ID Software is an automated genotyping software for forensic casework, databasing, and paternity data analysis.

After electrophoresis, the data collection software stores information for each sample in a .fsa file. Using GeneMapper™ ID Software v3.2.1 software, you can then analyze and interpret the data from the .fsa files.

#### Instruments

Refer to [“Instrument and software overview”](#) on page 16 for a list of compatible instruments.

**Before you start**

When using GeneMapper™ ID Software v3.2.1 to perform human identification (HID) analysis with AmpFℓSTR™ kits, be aware that:

- HID analysis requires at least one allelic ladder sample per run folder. Your laboratory can use multiple ladder samples in an analysis, provided individual laboratories conduct the appropriate validation studies.  
For multiple ladder samples, the GeneMapper™ ID Software calculates allelic bin offsets by using an average of all ladders that use the same panel within a run folder.
- Allelic ladder samples in an individual run folder are considered to be from a single run.  
When the software imports multiple run folders into a project, only the ladder(s) within their respective run folders are used for calculating allelic bin offsets and subsequent genotyping.
- Allelic ladder samples must be labeled as “Allelic Ladder” in the Sample Type column in a project. Failure to apply this setting for ladder samples results in failed analysis.
- Injections containing the allelic ladder must be analyzed with the same analysis method and parameter values that are used for samples to ensure proper allele calling.
- Alleles that are not in the AmpFℓSTR™ Allelic Ladders do exist. Off-ladder (OL) alleles may contain full and/or partial repeat units. An off-ladder allele is an allele that occurs outside the  $\pm 0.5$ -nt bin window of any known allelic ladder allele or virtual bin.  
**Note:** If a sample allele peak is called as an off-ladder allele, the sample result needs to be verified according to the laboratory’s protocol.

**Set up GeneMapper™ ID Software for data analysis****File names**

The file names shown in this section may differ from the file names you see when you download or import files. If you need help determining the correct files to use, contact your local Life Technologies Human Identification representative, or go to [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) ▶ **Software, Patches & Updates** ▶ **GeneMapper™ ID Software**.

**Workflow**

Before you can analyze sample (.fsa) files using GeneMapper™ ID Software v3.2.1 for the first time, you need to:

- Import panels and bins into the Panel Manager, as explained in “[Import panels and bins](#)” on page 35.
- Create an analysis method, as explained in , “[Create an analysis method](#)” on page 38.
- Create a size standard, as explained in “[Create size standard](#)” on page 43.
- Define custom views of analysis tables.  
Refer to Chapter 1 of the *GeneMapper™ ID Software Versions 3.1 and 3.2 Human Identification Analysis Tutorial* (Pub. no. 4335523) for more information.
- Define custom views of plots.  
Refer to Chapter 1 of the *GeneMapper™ ID Software Versions 3.1 and 3.2 Human Identification Analysis Tutorial* (Pub. no. 4335523) for more information.

## Import panels and bins

To import the SGM Plus™ panel and bin set into the GeneMapper™ ID Software v3.2.1 database:

1. Start the GeneMapper™ ID Software, then log in with the appropriate user name and password.

---

**IMPORTANT!** If you need logon instructions, refer to page 2-7 of the *GeneMapper™ ID Software Version 3.1 Human Identification Analysis User Guide* (Pub. no. 4338775).

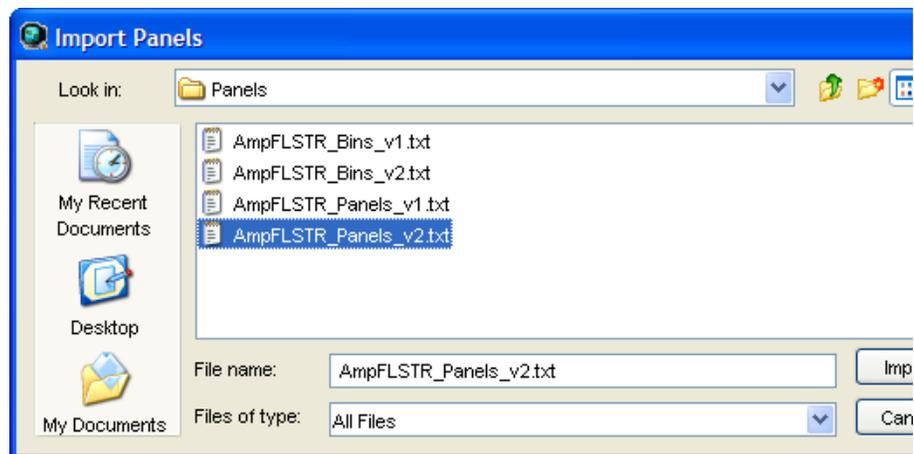
---

2. Select **Tools ▶ Panel Manager**.
3. Find, then open the folder containing the panels and bins:
  - a. Select **Panel Manager** in the navigation pane.



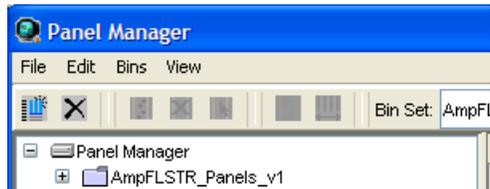
- b. Select **File ▶ Import Panels** to open the Import Panels dialog box.
  - c. Navigate to, then open the *x:\Applied Biosystems\GeneMapper\Panels* folder, where *x* is the drive on which the GeneMapper™ ID Software is installed.
4. Select **AmpFLSTR\_Panels\_v2.txt**, then click **Import**.

**Note:** Importing this file creates a new folder in the navigation pane of the Panel Manager, *AmpFLSTR\_Panels\_v2*. This folder contains the panels and associated markers.



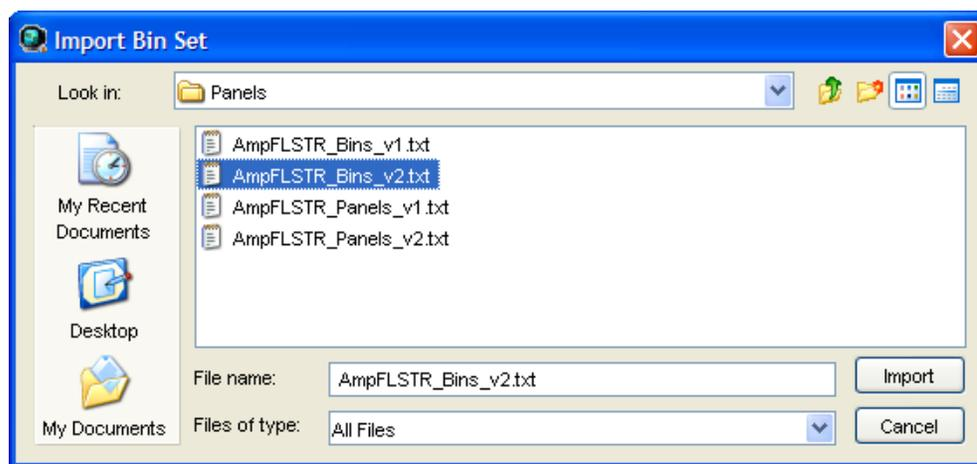
## 5. Import AmpFLSTR\_Bins\_v2.txt:

- a. Select the **AmpFLSTR\_Panels\_v2** folder in the navigation pane.



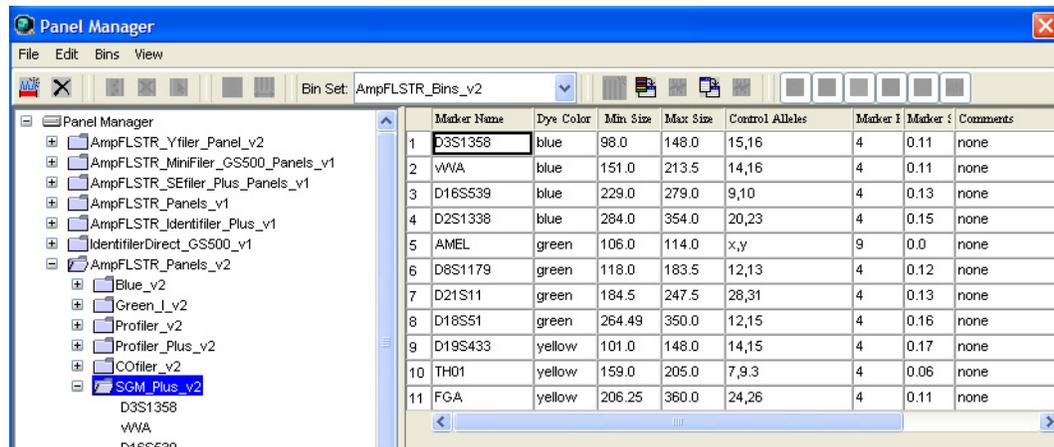
- b. Select **File** ► **Import Bin Set** to open the Import Bin Set dialog box.
- c. Navigate to, then open the *x:\Applied Biosystems\GeneMapper\Panels* folder.
- d. Select **AmpFLSTR\_Bins\_v2.txt**, then click **Import**.

**Note:** Importing this file associates the bin set with the panels in the AmpFLSTR\_Panels\_v2 folder.

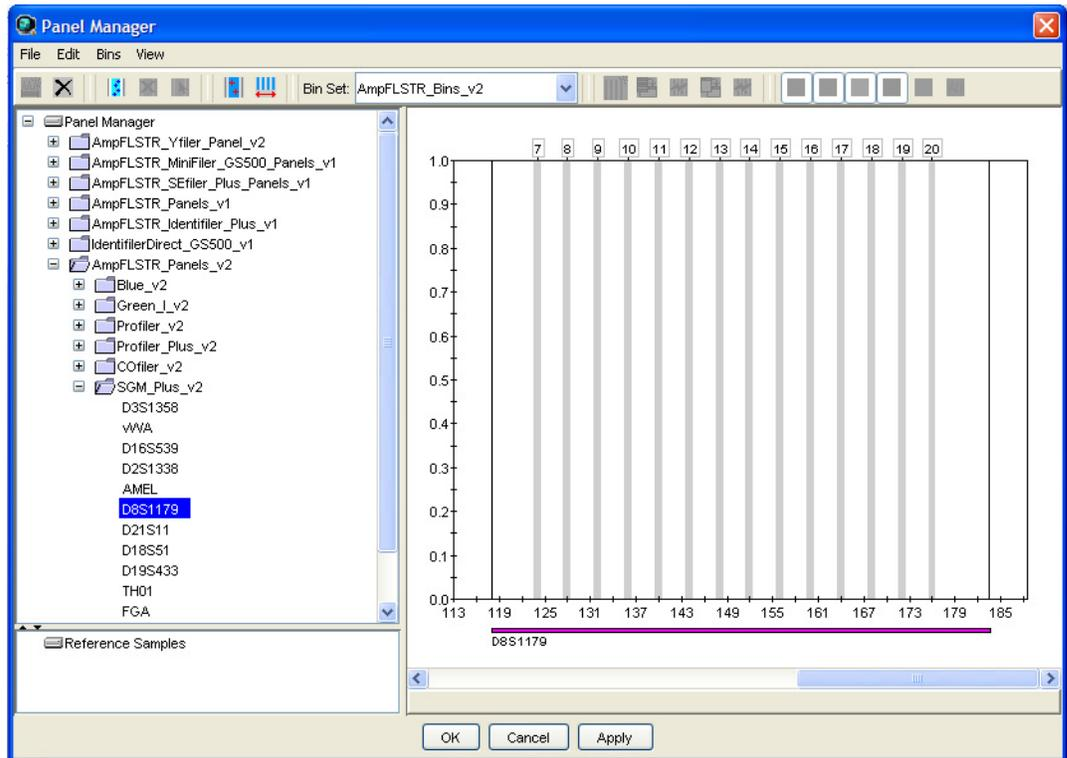


## 6. View the imported panels in the navigation pane:

- a. Double-click the **AmpFLSTR\_Panels\_v2** folder.
- b. Double-click the **SGM\_Plus\_v2** folder to display the panel information in the right pane and the markers below it.



7. View the markers and display the Bin view in the navigation pane:
  - a. Double-click the **AmpFLSTR\_Panels\_v2** folder to display its list of kits in the right pane.
  - b. Double-click the **SGM\_Plus\_v2** folder to display its list of markers below it.
  - c. Select **D8S1179** to display the Bin view for the marker in the right pane.



8. Click **Apply**, then **OK** to add the SGM\_Plus\_v2 panel and bin set to the GeneMapper™ ID Software database.

---

**IMPORTANT!** If you close the Panel Manager without clicking OK, the panels and bins are not imported into the GeneMapper™ ID Software database.

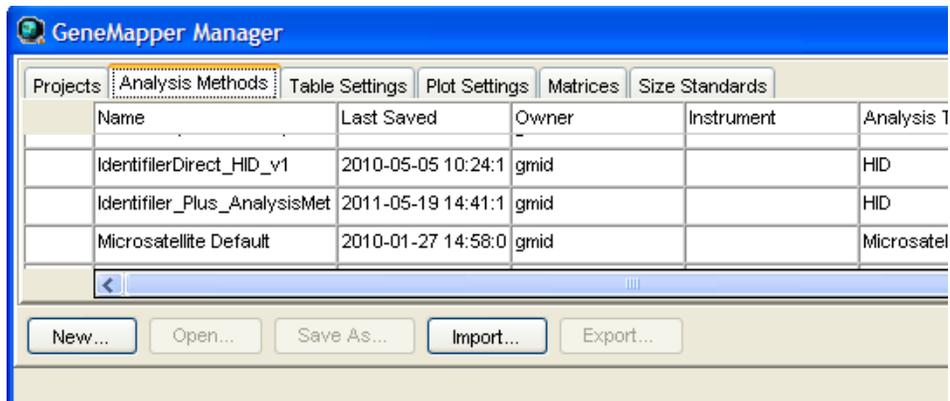
---

## Create an analysis method

The HID Advanced analysis method for the SGM Plus™ Kit uses the AmpFLSTR\_Bins\_v2 file described in [step 5 on page 36](#).

Use the following procedure to create a HID analysis method for the SGM Plus™ Kit.

1. Select **Tools** ▶ **GeneMapper Manager** to open the GeneMapper Manager.



2. Select the **Analysis Methods** tab, then click **New** to open the New Analysis Method dialog box.
3. Select **HID** and click **OK** to open the Analysis Method Editor with the General Tab selected.
4. The figures below show the settings for each tab of the Analysis Method Editor. Configure settings as shown unless the instructions state otherwise.
 

**Note:** The Analysis Method Editor closes when you save your settings. To complete this step quickly, do not save the analysis method until you finish entering settings in all of the tabs.
5. After you enter settings in all tabs, click **Save**.

## General tab settings

The screenshot shows the 'Analysis Method Editor - HID' dialog box with the 'General' tab selected. The 'Analysis Method Description' section contains the following fields:

- Name:** SGMPlus\_AnalysisMethod\_v1
- Description:** (Empty text area with scroll bars)
- Instrument:** (Empty text field)
- Analysis Type:** HID

At the bottom of the dialog are 'OK' and 'Cancel' buttons.

In the Name field, either type the name as shown, or enter a name of your choosing. The Description and Instrument fields are optional.

## Allele tab settings

Analysis Method Editor - HID

General | **Allele** | Peak Detector | Peak Quality | Quality Flags

Bin Set: AmpFLSTR\_Bins\_v2

Use marker-specific stutter ratio if available

Marker Repeat Type :		Tri	Tetra	Penta	Hexa
Cut-off Value		0.0	0.0	0.0	0.0
MinusA Ratio		0.0	0.0	0.0	0.0
MinusA Distance	From	0.0	0.0	0.0	0.0
	To	0.0	0.0	0.0	0.0
Minus Stutter Ratio		0.0	0.0	0.0	0.0
Minus Stutter Distance	From	0.0	3.25	0.0	0.0
	To	0.0	4.75	0.0	0.0
Plus Stutter Ratio		0.0	0.0	0.0	0.0
Plus Stutter Distance	From	0.0	0.0	0.0	0.0
	To	0.0	0.0	0.0	0.0

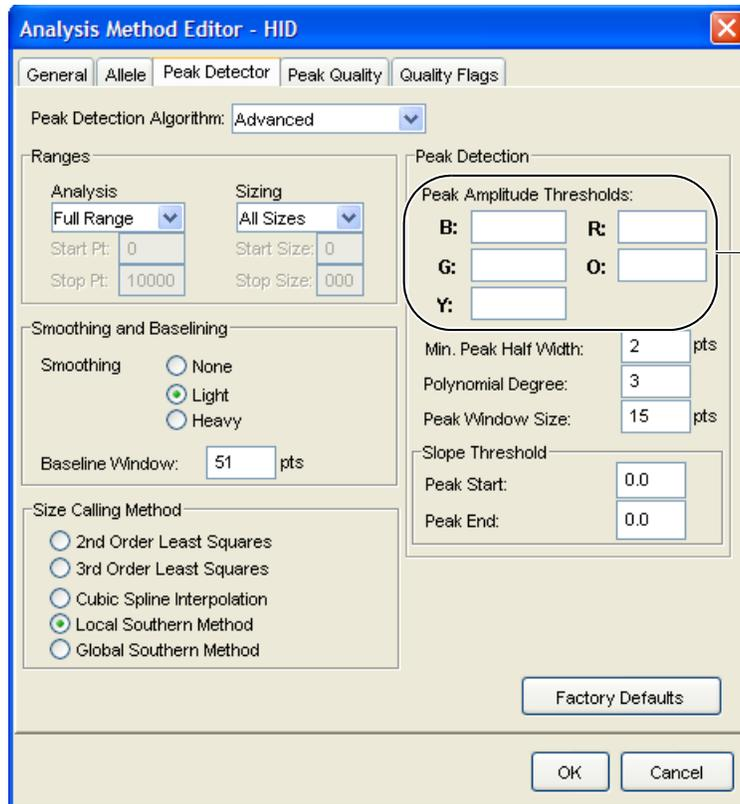
Amelogenin Cutoff: 0.0

Range Filter... Factory Defaults

OK Cancel

- In the Bin Set field, select the **AmpFLSTR\_Bins\_v2** bin set imported previously and configure the stutter distance parameters as shown.
- GeneMapper™ ID Software v3.2.1 allows you to specify four types of marker repeat motifs: tri, tetra, penta, and hexa. You can enter parameter values for each type of repeat in the appropriate column.
- Specify the stutter ratio:
  - To apply the stutter ratios listed in the Allele tab for single-source data, deselect the “Use marker-specific stutter ratio if available” check box (selected by default). Perform appropriate internal validation studies to determine the appropriate filter setting to use.  
**Note:** Applying global stutter ratios may reduce the editing required for single-source sample data.
  - To apply the stutter ratios contained in the AmpFLSTR\_Panels\_v2.txt file, select the “Use marker-specific stutter ratio if available” check box (selected by default). Perform appropriate internal validation studies to determine the appropriate filter setting to use.

Peak Detector tab settings



Perform internal validation studies to determine settings

**IMPORTANT!** Perform the appropriate internal validation studies to determine the peak amplitude thresholds for interpretation of SGM Plus™ Kit data.

Fields include:

- **Peak amplitude thresholds** – The software uses these parameters to specify the minimum peak height, in order to limit the number of detected peaks. Although GeneMapper™ ID Software displays peaks that fall below the specified amplitude in electropherograms, the software does not label or determine the genotype of these peaks.
- **Size calling method** – The SGM Plus™ Kit has been validated using the Local Southern sizing method. Before using other sizing methods, perform internal validation studies.

## Peak Quality tab settings

The screenshot shows the 'Analysis Method Editor - HID' window with the 'Peak Quality' tab selected. The 'Signal level' section is circled in red, containing three input fields: 'Homozygous min peak height', 'Heterozygous min peak height', and 'Min peak height ratio'. The 'Peak morphology' section has 'Max peak width (basepairs)' set to 1.5. The 'Pull-up peak' section has 'Pull-up ratio' set to 0.05. The 'Allele number' section has 'Max expected alleles' set to 2. At the bottom right, there are buttons for 'Factory Defaults', 'OK', and 'Cancel'.

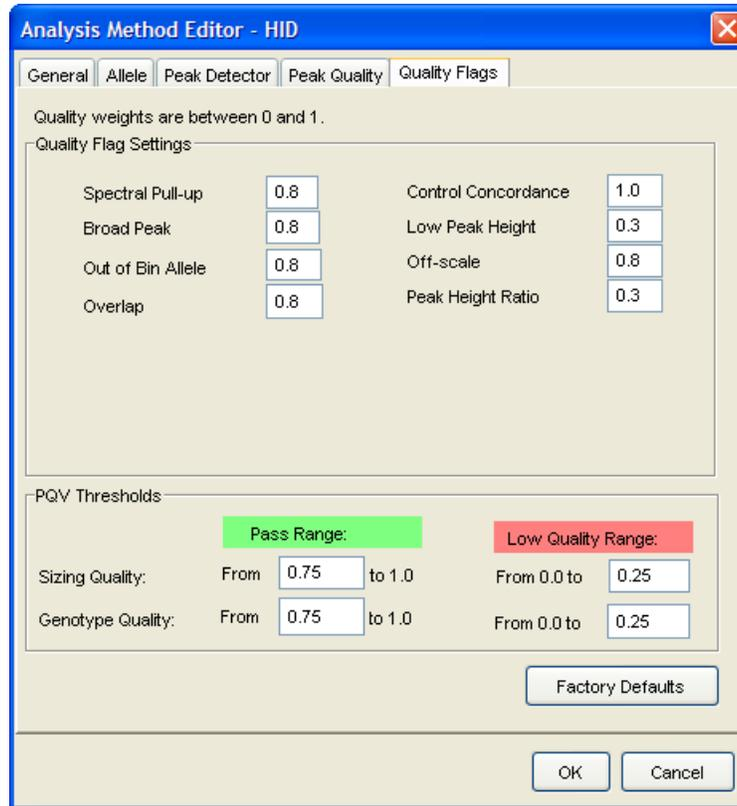
Perform internal validation studies to determine settings

---

**IMPORTANT!** Perform the appropriate internal validation studies to determine the minimum heterozygous and homozygous minimum peak height thresholds and the minimum peak height ratio threshold for interpretation of SGM Plus™ Kit data.

---

## Quality Flags tab settings



**IMPORTANT!** The values shown are the software defaults and are the values we used during developmental validation. Perform the appropriate internal validation studies to determine the appropriate values for interpretation of SGM Plus™ Kit data.

## Create size standard

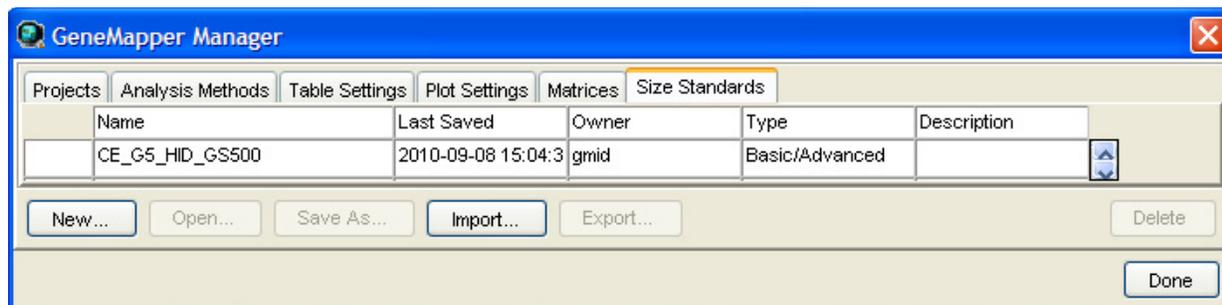
The GeneScan™ 500 ROX™ Size Standard for the SGM Plus™ Kit uses the following size standard peaks in its definitions: 75, 100, 139, 150, 160, 200, 300, 340, 350, 400, and 450.

**Note:** The 250-nt peak in the GeneScan™ 500 ROX™ Size Standard is not included in the size standard definition. This peak can be used as an indicator of precision within a run.

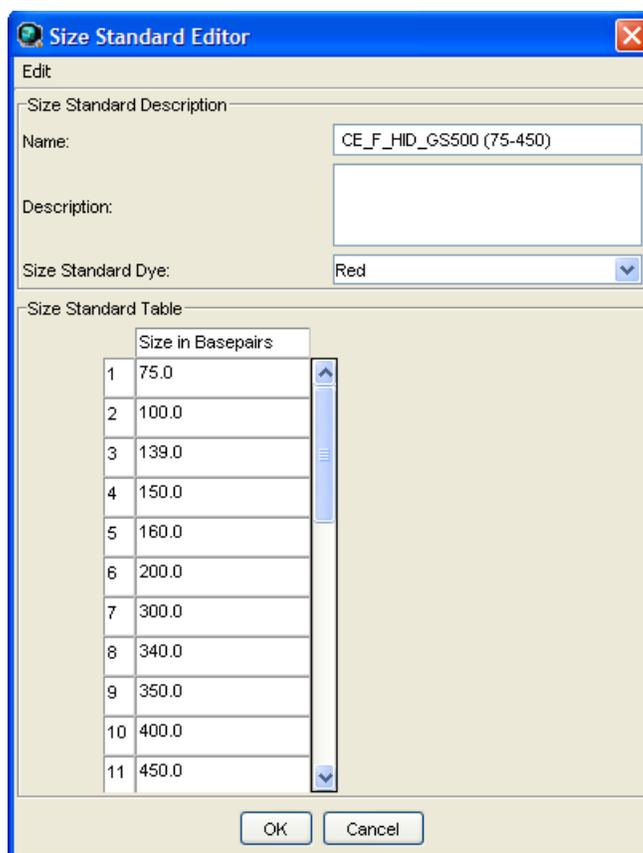
Use the following procedure to create the appropriate size standard:

1. Select **Tools ▶ GeneMapper Manager** to open the GeneMapper Manager.

2. Select the **Size Standards** tab, click **New**, select the **Basic or Advanced** radio button, then click **OK**.



3. Enter a name (for example, CE\_F\_HID\_GS500(75-450)). In the Size Standard Dye field, select **Red**. In the Size Standard Table, enter the sizes specified in [on page 43](#). The example below is for the GeneScan™ 500 ROX™ Size Standard.



## Analyze and edit sample files with GeneMapper™ ID Software

1. In the Project window, select **File ▶ Add Samples** to Project, then navigate to the disk or directory containing the sample files.
2. Apply analysis settings to the samples in the project.

Parameter	Settings
Sample Type	Select the sample type.
Analysis Method	SGMPlus_AnalysisMethod_v1 (or the name of the analysis method you created)
Panel	SGM_Plus_v2
Size Standard	CE_F_HID_GS500(75-450) <sup>†</sup> (or the name of the size standard you created)

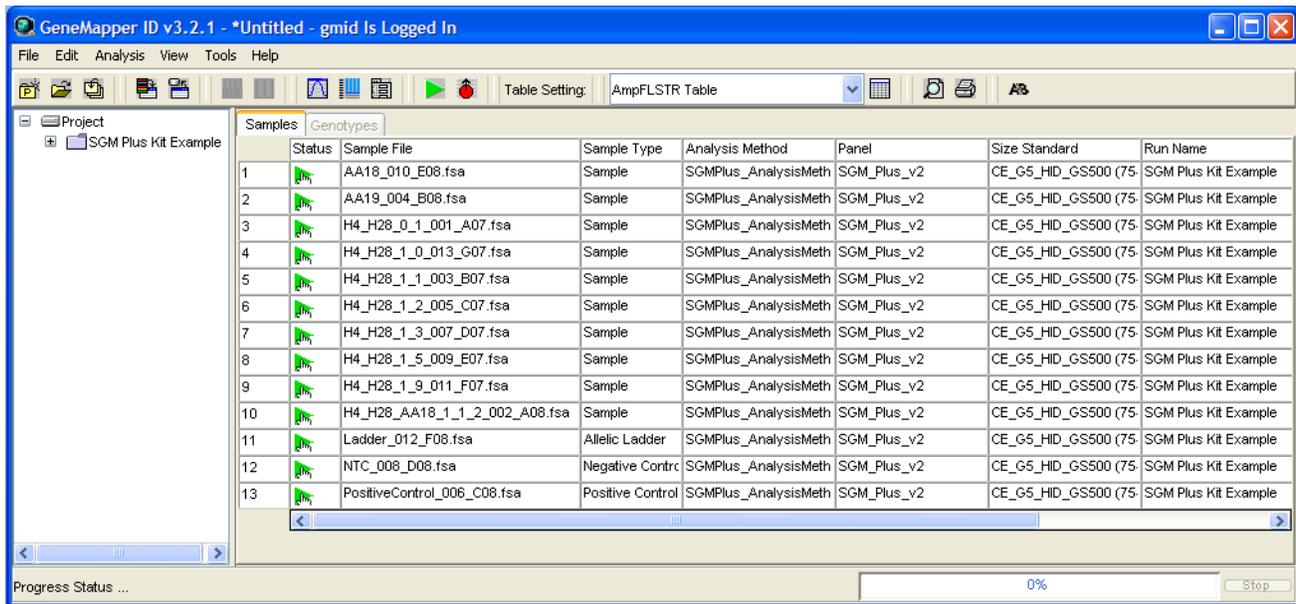
<sup>†</sup> The SGM Plus™ Kit was originally validated using the GeneScan™ 500 ROX™ Size Standard. If you use the GeneScan™ 400 HD Size Standard as an alternative, perform the appropriate internal validation studies to support the use of this size standard with the SGM Plus™ Kit.

**Note:** For more information about how the Size Caller works, refer to the *GeneScan™ Analysis Software for the Windows™ NT Operating System Overview of the Analysis Parameters and Size Caller User Bulletin* (Pub. no. 4335617).

3. Click  (Analyze), enter a name for the project (in the Save Project dialog box), then click **OK** to start analysis.
  - The status bar displays the progress of analysis:
    - As a completion bar extending to the right with the percentage indicated
    - With text messages on the left

- The table displays the row of the sample currently being analyzed in green (or red if analysis failed for the sample).
- The Genotypes tab becomes available after analysis (see the figure on the next page).

### Project window after analysis



For more information about any of these tasks, refer to the GeneMapper™ ID Software Version 3.1 Human Identification Analysis User Guide (Pub. no. 4338775).

## Examine and edit a project

You can display electropherogram plots from the Samples and Genotypes tabs of the Project window to examine the data. These procedures start with the Samples tab of the Project window (assuming the analysis is complete).

## For more information

For details about GeneMapper™ ID Software features, allele filters, peak detection algorithms, and project editing, refer to:

- *GeneMapper™ ID Software Versions 3.1 and 3.2 Human Identification Analysis Tutorial* (Pub. no. 4335523)
- *GeneMapper™ ID Software Version 3.1 Human Identification Analysis User Guide* (Pub. no. 4338775)
- *Installation Procedures and New Features for GeneMapper™ ID Software Software Version v3.2 User Bulletin* (Pub. no. 4352543)



## Section 4.2 GeneMapper™ ID-X Software

### Overview of GeneMapper™ ID-X Software

GeneMapper™ ID-X Software is an automated genotyping software for forensic casework, databasing, and paternity data analysis.

After electrophoresis, the data collection software stores information for each sample in a .fsa or .hid file. Using GeneMapper™ ID-X Software v1.0.1 or higher you can then analyze and interpret the data from the .fsa or .hid files.

#### Instruments

Refer to [“Instrument and software overview”](#) on page 16 for a list of compatible instruments.

#### Before you start

When using GeneMapper™ ID-X Software v1.0.1 or higher to perform human identification (HID) analysis with AmpF $\mathcal{L}$ STR™ kits, be aware that:

- HID analysis requires at least one allelic ladder sample per run folder. Your laboratory can use multiple ladder samples in an analysis, provided individual laboratories conduct the appropriate validation studies.  
For multiple ladder samples, the GeneMapper™ ID-X Software calculates allelic bin offsets by using an average of all ladders that use the same panel within a run folder.
- Allelic ladder samples in an individual run folder are considered to be from a single run.  
When the software imports multiple run folders into a project, only the ladder(s) within their respective run folders are used for calculating allelic bin offsets and subsequent genotyping.
- Allelic ladder samples must be labeled as “Allelic Ladder” in the Sample Type column in a project. Failure to apply this setting for ladder samples results in failed analysis.
- Injections containing the allelic ladder must be analyzed with the same analysis method and parameter values that are used for samples to ensure proper allele calling.
- Alleles that are not in the AmpF $\mathcal{L}$ STR™ Allelic Ladders do exist. Off-ladder (OL) alleles may contain full and/or partial repeat units. An off-ladder allele is an allele that occurs outside the  $\pm 0.5$ -nt bin window of any known allelic ladder allele or virtual bin.  
**Note:** If a sample allele peak is called as an off-ladder allele, the sample result needs to be verified according to the laboratory’s protocol.

## Set up GeneMapper™ ID-X Software for data analysis

### Panel, bin, and stutter file version

The file names shown in this section may differ from the file names you see when you download or import files. If you need help determining the correct files to use, contact your local Life Technologies Human Identification representative, or go to [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) ▶ **Software, Patches & Updates** ▶ **GeneMapper™ ID-X Software**.

The instructions and examples in this section refer to the latest version of panel, bin, and stutter file available at the time of publication.

### Before using the software for the first time

Before you use GeneMapper™ ID-X Software (v1.0.1 or higher for .fsa files, v1.2 or higher for .hid files) to analyze data for the first time, you must do the following:

1. Check the version of panel, bin, and stutter files installed with the GeneMapper™ ID-X Software as explained in “[Check panel, bin, and stutter file version](#)” below.
2. Check [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) ▶ **Software, Patches & Updates** ▶ **GeneMapper™ ID-X Software** to determine if newer files are available.
3. If updated files are available, download and import the files into the GeneMapper™ ID-X Software, as explained in “[Import panels, bins, and marker stutter](#)” on page 50.  
**Note:** When downloading new versions of analysis files, refer to the associated Read Me file for details of changes between software file versions. If you have validated previous file versions for data analysis, conduct the appropriate internal verification studies before using new file versions for operational analysis.
4. Create an analysis method, as explained in “[Create an analysis method](#)” on page 55.
5. Define custom views of analysis tables.  
Refer to Chapter 1 of the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Pub. no. 4375574) for more information.
6. Define custom views of plots.  
Refer to Chapter 1 of the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Pub. no. 4375574) for more information.

### Check panel, bin, and stutter file version

1. Start the GeneMapper™ ID-X Software, then log in with the appropriate user name and password.

---

**IMPORTANT!** For logon instructions, refer to the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Pub. no. 4375574).

---

2. Select **Tools** ▶ **Panel Manager**.

3. Check the version of files imported into the Panel Manager:

- a. Select **Panel Manager** in the navigation pane.
- b. Expand the Panel Manager folder and any sub-folders to identify the analysis file version already installed for your kit choice.



4. Check the version of files available for import into the Panel Manager:

- a. Select **Panel Manager**, then select **File ▶ Import Panels** to open the Import Panels dialog box.
  - b. Navigate to, then open the Panels folder and check the version of panel, bin, and stutter files installed.
5. If newer versions are available on the website, download and import as described below.

### Import panels, bins, and marker stutter

To import the SGM Plus™ Kit panel, bin set, and marker stutter from our web site into the GeneMapper™ ID-X Software database:

1. Download and open the file containing panels, bins, and marker stutter:
  - a. Go to [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) ▶ **Software, Patches & Updates ▶ GeneMapper™ ID-X Software**. Download the file **AmpFLSTR Analysis Files GMIDX**.
  - b. Unzip the file.
2. Start the GeneMapper™ ID-X Software, then log in with the appropriate user name and password.

---

**IMPORTANT!** For logon instructions, refer to the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Pub. no. 4375574).

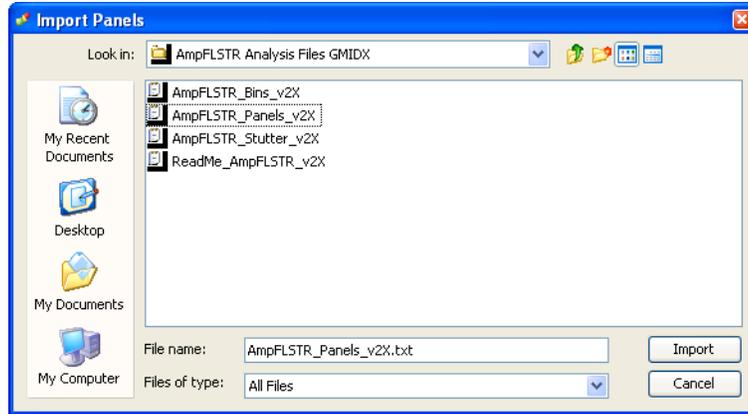
---

3. Select **Tools ▶ Panel Manager**.
4. Find, then open the folder containing the panels, bins, and marker stutter:
  - a. Select **Panel Manager** in the navigation pane.
  - b. Select **File ▶ Import Panels** to open the Import Panels dialog box.
  - c. Navigate to, then open the **AmpFLSTR Analysis Files GMIDX** folder that you unzipped in [step 1 on page 50](#).

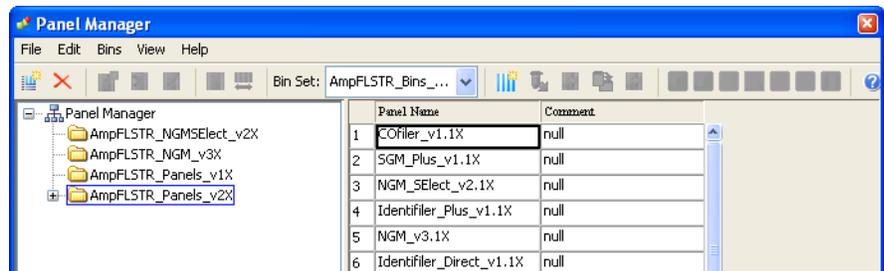


5. Select **AmpFLSTR\_Panels\_v2X** (or the version you installed), then click **Import**.

**Note:** Importing this file creates a new folder in the navigation pane of the Panel Manager “AmpFLSTR\_Panels\_v2X”. This folder contains panels for multiple AmpFLSTR™ kits and associated markers.



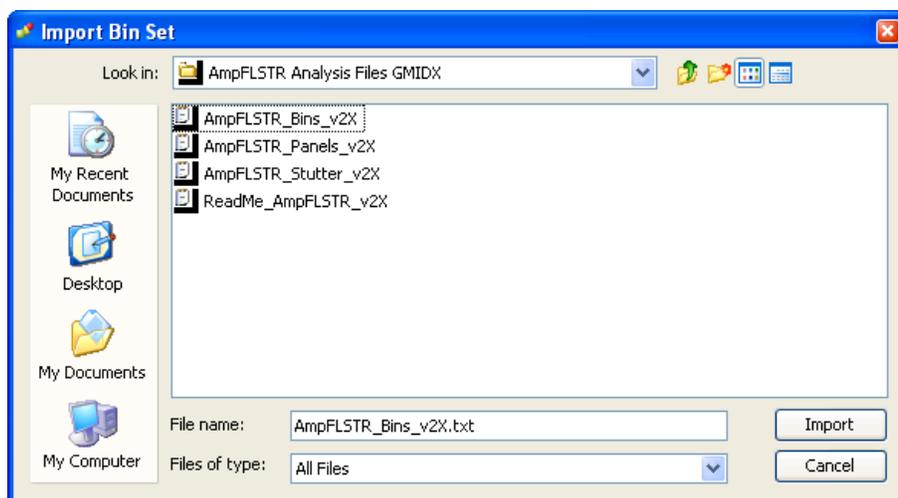
6. Import AmpFLSTR\_Bins\_v2X.txt:
  - a. Select the **AmpFLSTR\_Panels\_v2X** folder in the navigation pane.



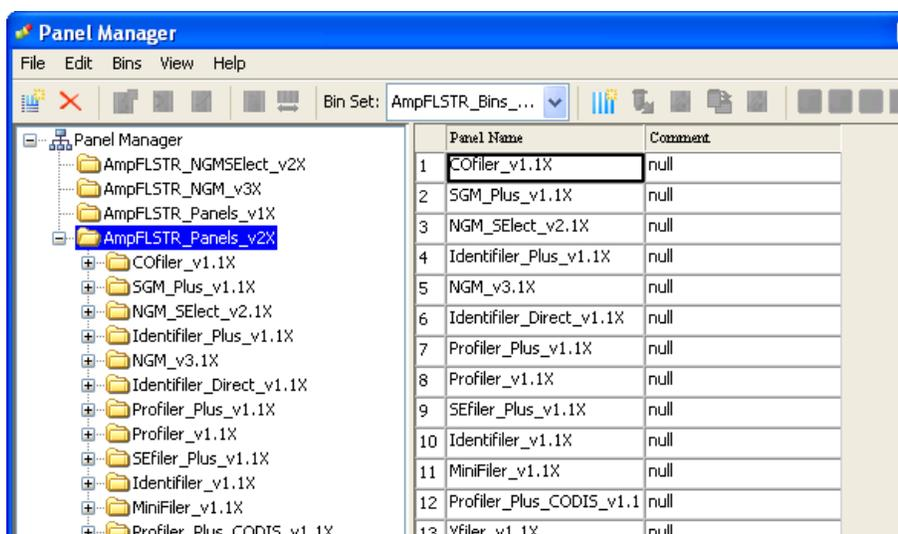
- b. Select **File ► Import Bin Set** to open the Import Bin Set dialog box.
  - c. Navigate to, then open the **AmpFLSTR Analysis Files GMIDX** folder.

- d. Select **AmpFLSTR\_Bins\_v2X.txt**, then click **Import**.

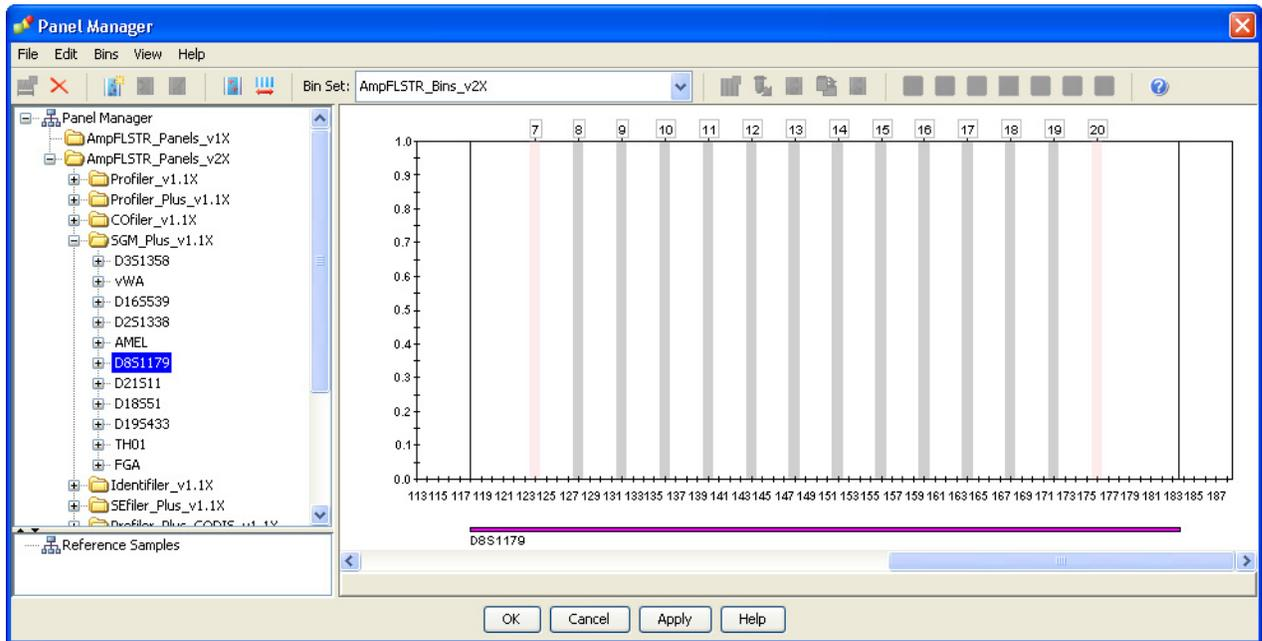
**Note:** Importing this file associates the bin set with the panels in the AmpFLSTR\_Panels\_v2X folder.



7. View the imported panels in the navigation pane:
- Double-click the **AmpFLSTR\_Panels\_v2X** folder.
  - Double-click the **SGM\_Plus\_v1.1X** folder to display the panel information in the right pane.

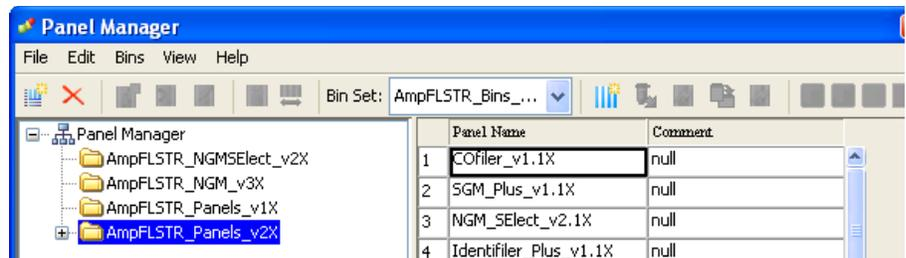


8. Select and expand **SGM\_Plus\_v1.1X** in the navigation pane, then select **D8S1179** to display the Bin view for the marker in the right pane.



9. Import AmpFLSTR\_Stutter\_v2X:

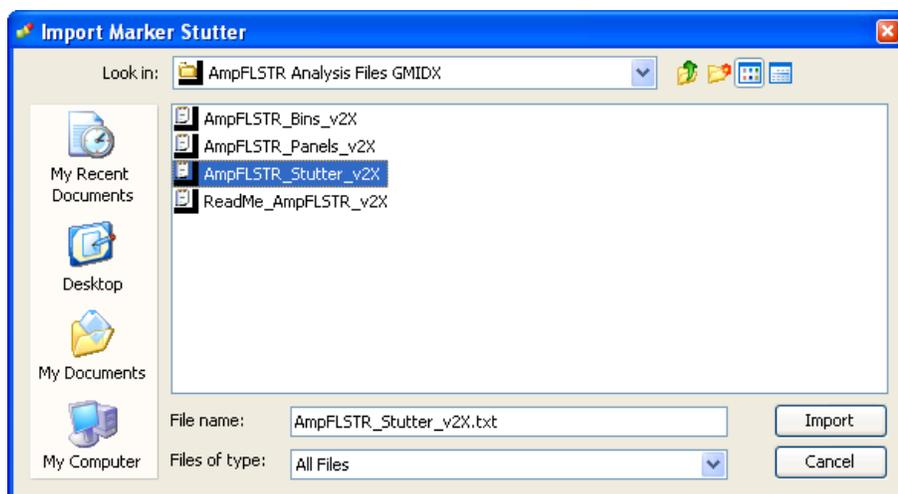
- a. Select the **AmpFLSTR\_Panels\_v2X** folder in the navigation panel.



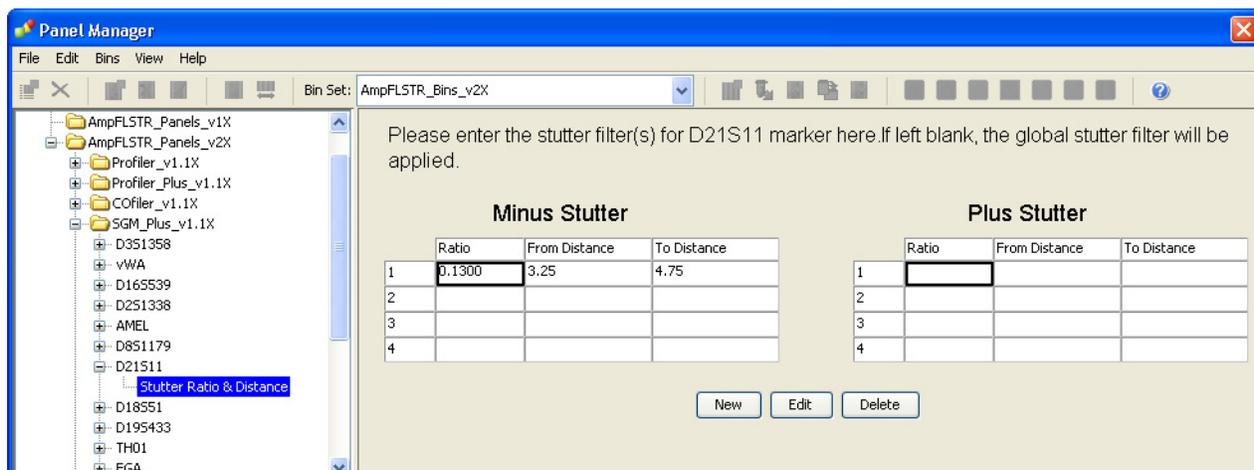
- b. Select **File** ► **Import Marker Stutter** to open the Import Marker Stutter dialog box.
- c. Navigate to, then open the **AmpFLSTR Analysis Files GMIDX** folder.

- d. Select **AmpFLSTR\_Stutter\_v2X**, then click **Import**.

**Note:** Importing this file associates the marker stutter ratio with the bin set in the AmpFLSTR\_Panels\_v2X folder.



10. View the imported marker stutters in the navigation pane:
- Double-click the **AmpFLSTR\_Panels\_v2X** folder to display its list of kits in the right pane.
  - Double-click the **SGM\_Plus\_v1.1X** folder to display its list of markers below it.
  - Double-click **D21S11** to display the Stutter Ratio & Distance view for the marker in the right pane.



11. Click **Apply**, then **OK** to add the SGM Plus™ Kit panel, bin set, and marker stutter to the GeneMapper™ ID-X Software database.

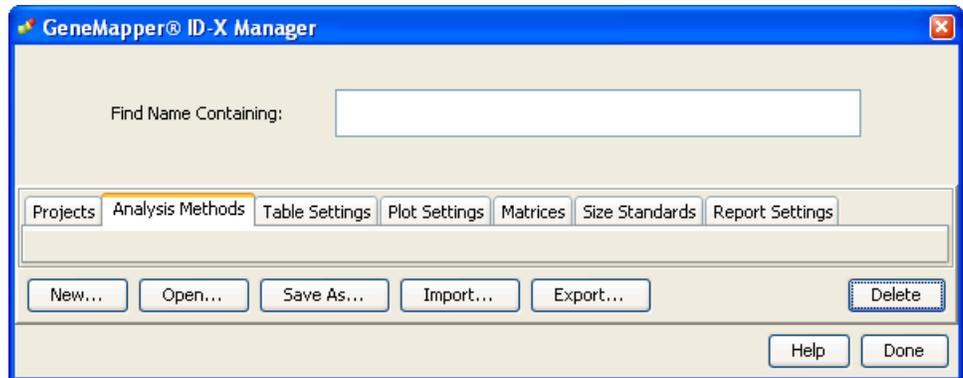
**IMPORTANT!** If you close the Panel Manager without clicking **Apply**, the panels, bin sets, and marker stutter will not be imported into the GeneMapper™ ID-X Software database.

## Create an analysis method

Use the following procedure to create an analysis method for the SGM Plus™ Kit.

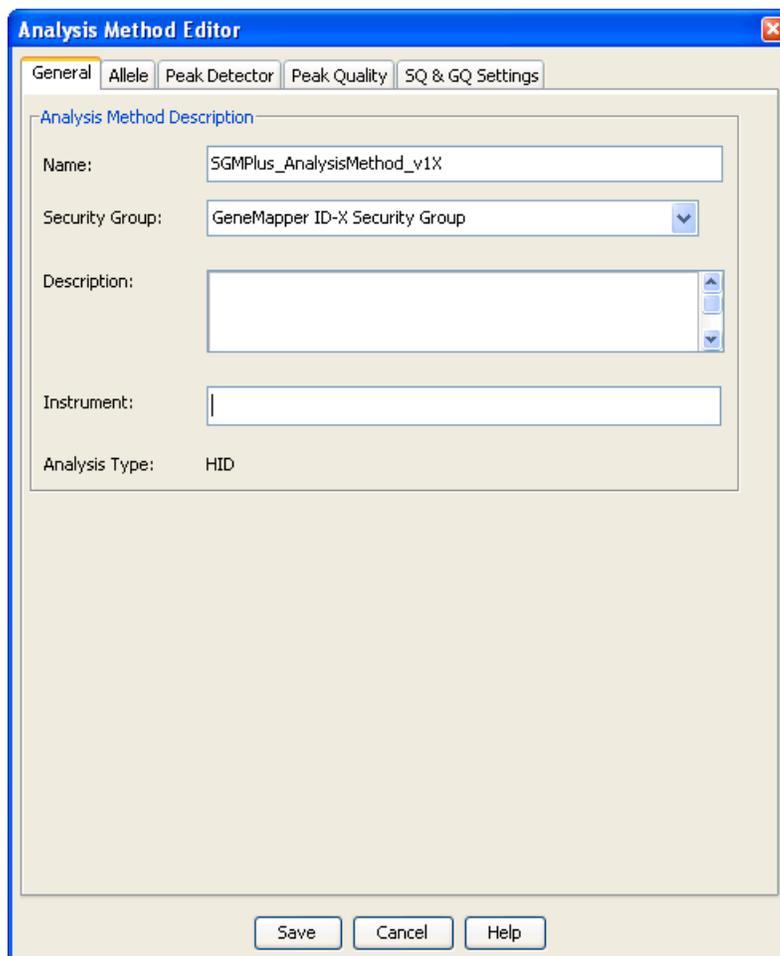
**IMPORTANT!** Analysis methods are version-specific, so you must create an analysis method for each version of the software. For example, an analysis method created for GeneMapper™ ID-X version 1.2 is not compatible with earlier versions of GeneMapper™ ID-X Software or with GeneMapper™ ID Software version 3.2.1.

1. Select **Tools** ▶ **GeneMapper™ ID-X Manager** to open the GeneMapper™ ID-X Manager.



2. Select the **Analysis Methods** tab, then click **New** to open the Analysis Method Editor with the **General** tab selected.  
The figures below show the settings for each tab of the Analysis Method Editor. Configure the Analysis Method Editor tab settings as shown in the figures below, unless the instructions state otherwise.  
**Note:** The Analysis Method Editor closes when you save your settings. To complete this step quickly, do not save the analysis method until you finish entering settings in all of the tabs.
3. After you enter settings in all tabs, click **Save**.

## General tab settings



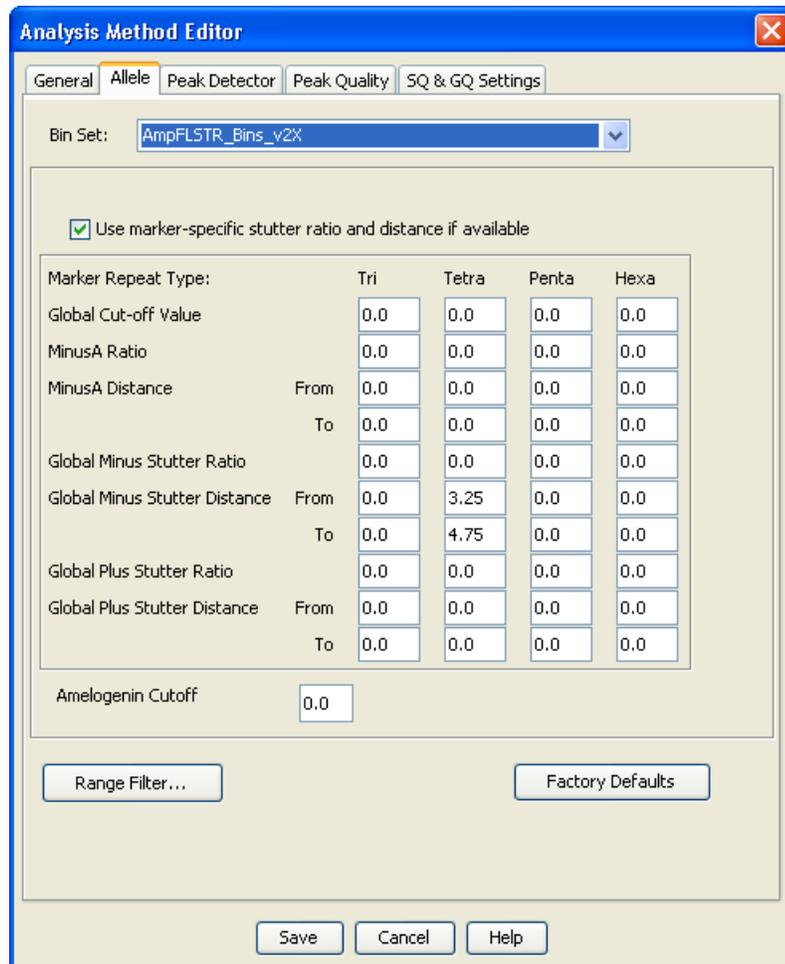
The screenshot shows the 'Analysis Method Editor' dialog box with the 'General' tab selected. The dialog has a title bar with a close button. Below the title bar are five tabs: 'General', 'Allele', 'Peak Detector', 'Peak Quality', and 'SQ & GQ Settings'. The 'General' tab is active and contains the following fields:

- Name:** A text box containing 'SGMPlus\_AnalysisMethod\_v1X'.
- Security Group:** A dropdown menu with 'GeneMapper ID-X Security Group' selected.
- Description:** A large text area with a vertical scrollbar, currently empty.
- Instrument:** A text box, currently empty.
- Analysis Type:** A label with the value 'HID'.

At the bottom of the dialog are three buttons: 'Save', 'Cancel', and 'Help'.

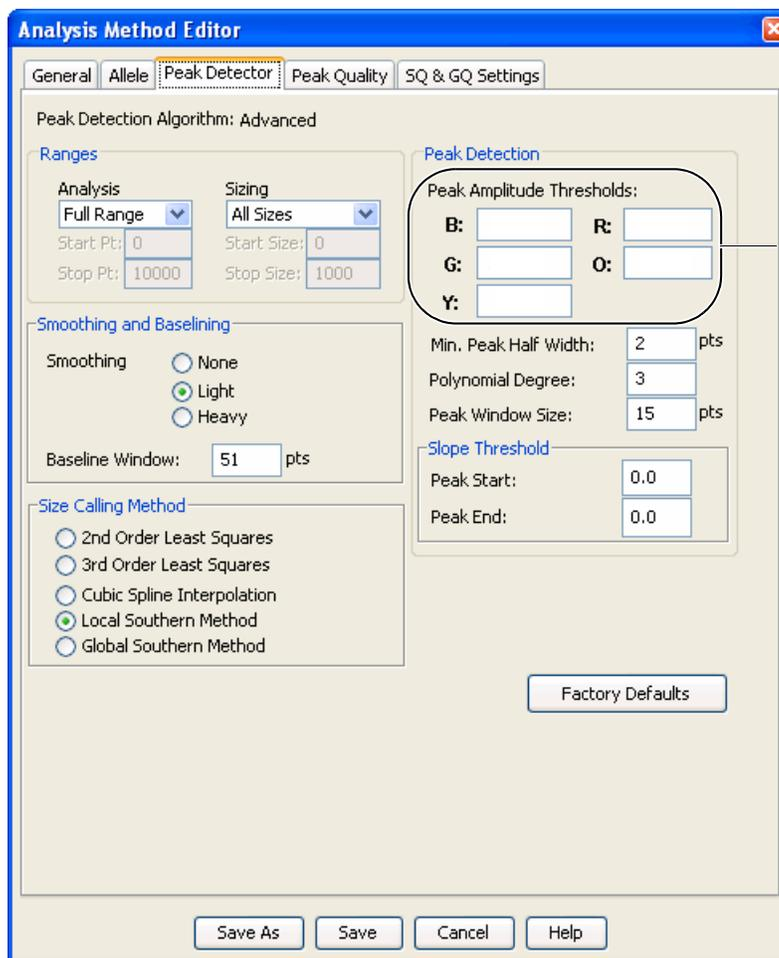
In the Name field, either type the name as shown or enter a name of your choosing. In the Security Group field, select the Security Group appropriate to your software configuration from the dropdown list. The Description and Instrument fields are optional.

## Allele tab settings



- In the Bin Set field, select the **AmpFLSTR\_Bins\_v2X** bin set and configure the stutter distance parameters as shown.
- GeneMapper™ ID-X Software v1.0.1 or higher allows you to specify 4 types of marker repeat motifs: tri, tetra, penta and hexa. You can enter parameter values for each type of repeat in the appropriate column.
- Specify the stutter ratio:
  - To apply the stutter ratios listed in the Allele tab for single-source data, deselect the “Use marker-specific stutter ratio if available” check box (selected by default). Perform appropriate internal validation studies to determine the appropriate filter setting to use.  
**Note:** Applying global stutter ratios may reduce the editing required for single-source sample data.
  - To apply the stutter ratios contained in the AmpFLSTR\_Stutter\_v2X file, select the “Use marker-specific stutter ratio if available” check box (selected by default). Perform appropriate internal validation studies to determine the appropriate filter setting to use.

## Peak Detector tab settings



Perform internal validation studies to determine settings

---

**IMPORTANT!** Perform the appropriate internal validation studies to determine the appropriate peak amplitude thresholds for interpretation of SGM Plus™ Kit data.

---

Fields include:

- **Peak amplitude thresholds** – The software uses these parameters to specify the minimum peak height, in order to limit the number of detected peaks. Although GeneMapper™ ID-X Software displays peaks that fall below the specified amplitude in electropherograms, the software does not label or determine the genotype of these peaks.
- **Size calling method** – The SGM Plus™ Kit has been validated using the Local Southern sizing method. Before using other sizing methods, perform internal validation studies.
- **Normalization** – A Normalization checkbox is available on this tab in GeneMapper™ ID-X Software v1.2 for use in conjunction with data run on the Applied Biosystems 3500 Series Genetic Analyzers. Normalization cannot be applied to 4-dye data, so this feature is not for use with SGM Plus™ Kit data.

## Peak Quality tab settings

Analysis Method Editor

General Allele Peak Detector **Peak Quality** SQ & GQ Settings

Min/Max Peak Height (LPH/MPH)

Homozygous min peak height

Heterozygous min peak height

Max Peak Height (MPH)

Peak Height Ratio (PHR)

Min peak height ratio

Broad Peak (BD)

Max peak width (basepairs)

Allele Number (AN)

Max expected alleles

Allelic Ladder Spike

Cut-off Value

Factory Defaults

Save Cancel Help

Perform internal validation studies to determine settings

---

**IMPORTANT!** Perform the appropriate internal validation studies to determine the minimum heterozygous and homozygous minimum peak height thresholds, maximum peak height threshold and the minimum peak height ratio threshold for interpretation of SGM Plus™ Kit data.

---

## SQ & GQ tab settings

Analysis Method Editor

General Allele Peak Detector Peak Quality **SQ & GQ Settings**

Quality weights are between 0 and 1.

Sample and Control GQ Weighting

Broad Peak (BD)	0.8	Allele Number (AN)	1.0
Out of Bin Allele (BIN)	0.8	Low Peak Height (LPH)	0.3
Overlap (OVL)	0.8	Max Peak Height (MPH)	0.3
Marker Spike (SPK)	0.3	Off-scale (OS)	0.8
		Peak Height Ratio (PHR)	0.3

Control Concordance (CC) Weight = 1.0 (Only applicable to controls)

SQ Weighting

Broad Peak (BD)	0.5
-----------------	-----

Allelic Ladder GQ Weighting

Spike (SSPK/SPK)	1	Off-scale (OS)	1
------------------	---	----------------	---

SQ & GQ Ranges

	<b>Pass Range:</b>	<b>Low Quality Range:</b>
Sizing Quality:	From 0.75 to 1.0	From 0.0 to 0.25
Genotype Quality:	From 0.75 to 1.0	From 0.0 to 0.25

Reset Defaults

Save As Save Cancel Help

**IMPORTANT!** The values shown are the software defaults and are the values we used during developmental validation. Perform appropriate internal validation studies to determine the appropriate values to use.

## Create size standard (optional)

The CE\_F\_HID\_GS500 (75–450) size standard definition is installed with the software for use with the SGM Plus™ Kit and contains the following size standard peaks:

### GeneScan™ 500 ROX™ Size Standard

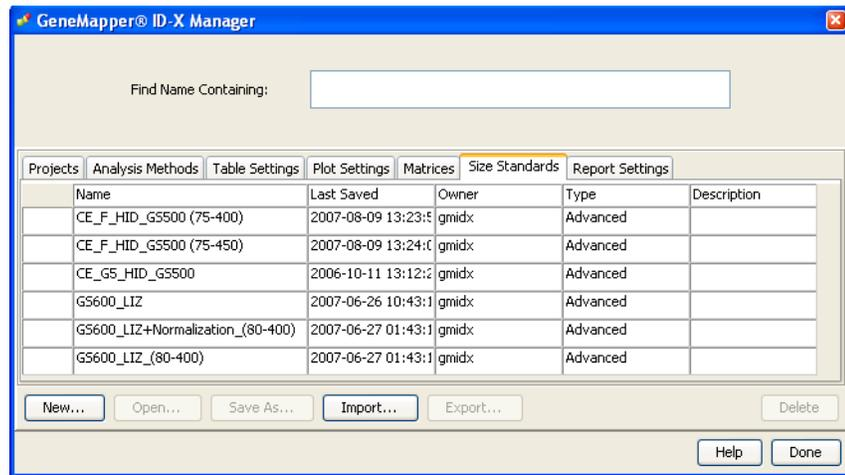
75, 100, 139, 150, 160, 200, 300, 340, 350, 400, and 450

**Note:** The 250-nt peak in the GeneScan™ 500 ROX™ Size Standard is not included in the size standard definition. This peak can be used as an indicator of precision within a run.

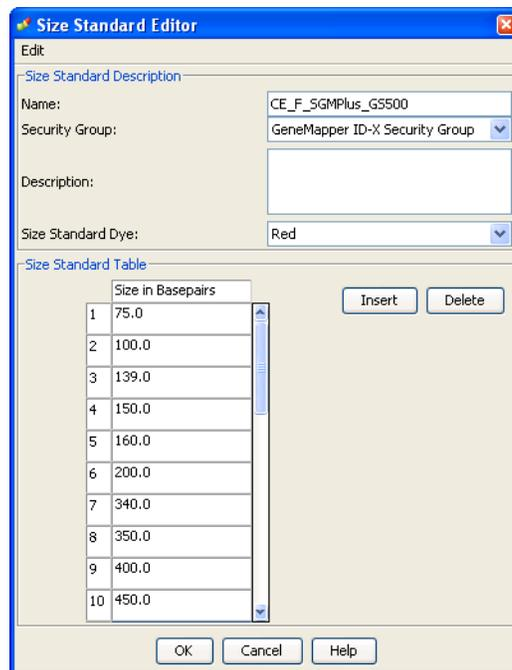
Use the following procedure if you want to create your own size standard:

1. Select **Tools** ▶ **GeneMapper Manager** to open the GeneMapper Manager.

2. Select the **Size Standards** tab, then click **New**.



3. Enter a name. In the Size Standard Dye field, select **Red**. In the Size Standard Table, enter the sizes specified in [on page 60](#). The example below is for the GeneScan™ 500 ROX™ Size Standard.



## Analyze and edit sample files with GeneMapper™ ID-X Software

1. In the Project window, select **File ▶ Add Samples to Project**, then navigate to the disk or directory containing the sample files.
2. Apply analysis settings to the samples in the project.

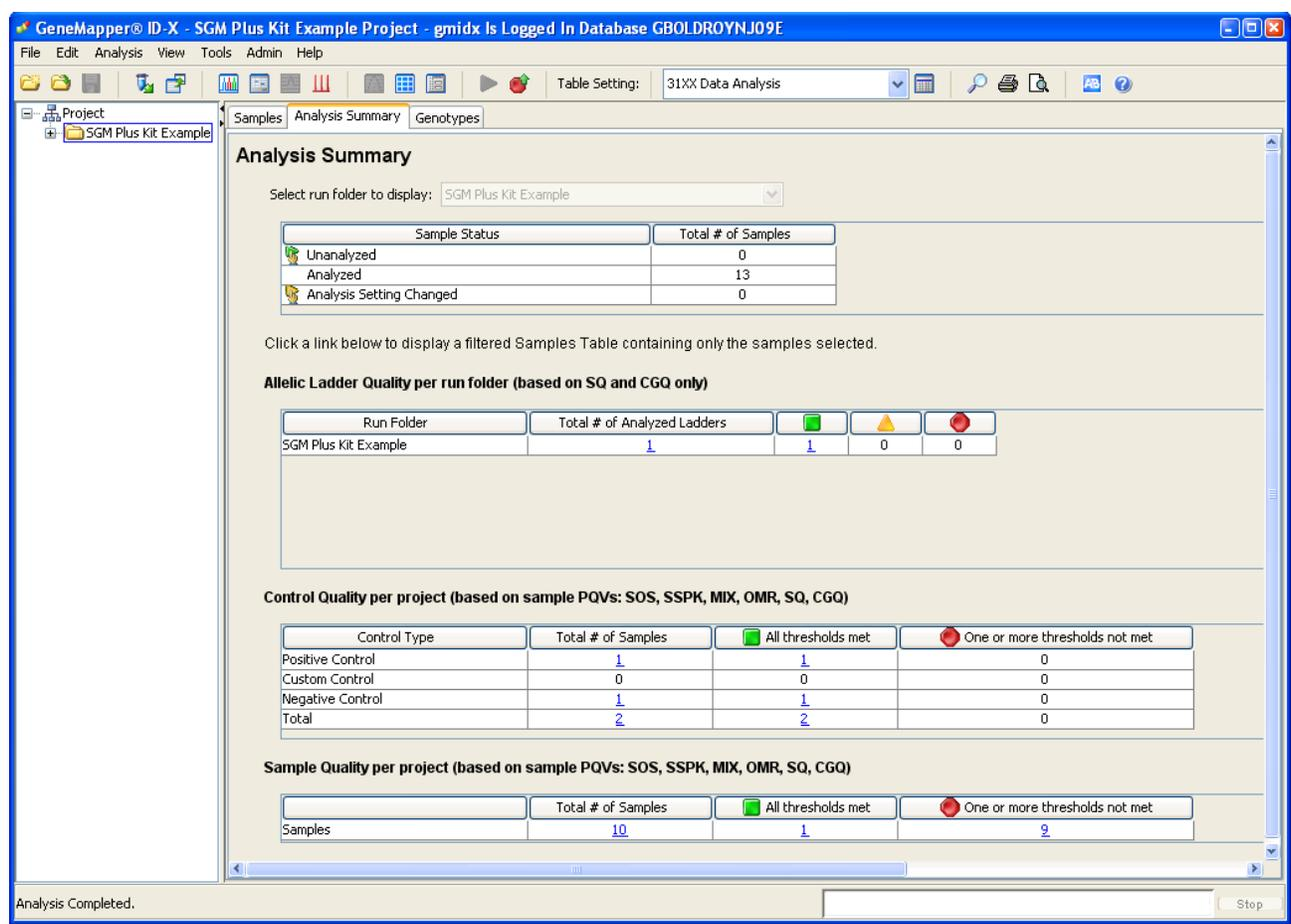
Parameter	Settings
Sample Type	Select the sample type.
Analysis Method	SGMPlus_AnalysisMethod_v1X (or the name of the analysis method you created)
Panel	SGM_Plus_v1.1X
Size Standard	CE_F_GS500(75-450) <sup>†</sup> (or the name of the size standard you created)

† The SGM Plus™ Kit was originally validated using the GeneScan™ 500 ROX™ Size Standard. If you use the GeneScan™ 400 HD Size Standard as an alternative, perform the appropriate internal validation studies to support the use of this size standard with the SGM Plus™ Kit.

**Note:** For more information about how the Size Caller works, refer to the *GeneScan™ Analysis Software for the Windows™ NT Operating System Overview of the Analysis Parameters and Size Caller User Bulletin* (Pub. no. 4335617).

3. Click  (**Analyze**), enter a name for the project (in the Save Project dialog box), then click **OK** to start analysis.
  - The status bar displays the progress of analysis as a completion bar extending to the right with the percentage indicated.
  - The table displays the row of the sample currently being analyzed in green (or red if analysis failed for the sample).
  - The Analysis Summary tab is displayed and the Genotypes tab becomes available upon completion of the analysis.

**Analysis summary window after analysis**



## Examine and edit a project

You can display electropherogram plots from the Samples and Genotypes tabs of the Project window to examine the data. These procedures start with the Analysis Summary tab of the Project window (assuming the analysis is complete).

## For more information

For more information, refer to:

- *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Pub. no. 4375574)
- *GeneMapper™ ID-X Software Version 1.0 Quick Reference Guide* (Pub. no. 4375670)
- *GeneMapper™ ID-X Software Version 1.0 Reference Guide* (Pub. no. 4375671)
- *GeneMapper™ ID-X Software Version 1.1(Mixture Analysis) Getting Started Guide* (Pub. no. 4396773)
- *GeneMapper™ ID-X Software Version 1.2 Reference Guide* (Pub. no. 4426481)
- *GeneMapper™ ID-X Software Version 1.2 Quick Reference Guide* (Pub. no. 4426482)

■ Section 5.1 Developmental Validation .....	66
Overview .....	66
Developmental validation.....	66
Accuracy, reproducibility, and precision.....	67
Extra peaks in the electropherogram.....	74
Characterization of loci .....	82
Species specificity .....	83
Sensitivity .....	85
Stability .....	86
Population data.....	92
Mutation rate.....	99
Probability of identity .....	100
Probability of paternity exclusion .....	100
■ Section 5.2 Performance Validation After Buffer and Enzyme Component Replacement .....	102
Overview .....	102
Experiments.....	102
Reproducibility study .....	103
Sensitivity study .....	106
Inhibition study .....	108
Conclusions.....	112

## Section 5.1 Developmental Validation

### Overview

This section provides results of the developmental validation experiments we performed using the SGM Plus™ Kit.

These studies meet and exceed those recommended in the Technical Working Group on DNA Analysis Methods (TWGDAM) guidelines as well as the DNA Advisory Board (DAB) Quality Assurance Standards, effective October 1, 1998 (Technical Working Group on DNA Analysis Methods, 1995; DNA Advisory Board, Federal Bureau of Investigation, U.S. Department of Justice, 1998). These studies also address the guidelines outlined in the ENFSI DNA Working Group Quality Assurance Programme for DNA Laboratories.

### Importance of validation

Validation of a DNA typing procedure for human identification applications is an evaluation of the procedure's efficiency, reliability, and performance characteristics. By challenging the procedure with samples commonly encountered in forensic and parentage laboratories, the validation process uncovers attributes and limitations which are critical for sound data interpretation in casework (Sparkes, Kimpton, Watson *et al.*, 1996; Sparkes, Kimpton, Gilbard *et al.*, 1996; Wallin *et al.*, 1998).

### Experiment conditions

This chapter discusses many of the experiments we performed and provides examples of results obtained. We chose conditions that produced optimum PCR product yield and that met reproducible performance standards. It is our opinion that while these experiments are not exhaustive, they are appropriate for a manufacturer of STR kits intended for forensic and/or parentage testing use. Each laboratory using the SGM Plus™ Kit should perform their own internal validation studies.

## Developmental validation

### DAB 8.1.1 Developmental Validation

*"Developmental validation that is conducted shall be appropriately documented."* (DNA Advisory Board, 1998).

Critical reagent concentrations and reaction conditions (such as thermal cycling parameters, AmpliTaq Gold™ DNA polymerase activation, cycle number) to produce reliable, locus-specific amplification and appropriate sensitivity have been determined.

### PCR components

The concentration of each component of the SGM Plus™ Kit was examined. The PCR components are Tris-HCl (pH 8.3), KCl, dNTPs, primers, AmpliTaq Gold™ DNA Polymerase, MgCl<sub>2</sub>, bovine serum albumin, and sodium azide. The concentration for a particular component was established to be in the window that meets the reproducible performance characteristics of specificity and sensitivity.

After the optimal concentration was determined for a single component, the others were tested sequentially until it was determined that each component was at the optimal concentration relative to the concentrations of the other components in the master mix. The optimized SGM Plus™ Kit provides the required degree of specificity such that it is specific to primates for the species tested (with the exception of the amelogenin locus, see “Nonhuman studies” on page 84) and does not produce nonspecific mispriming artifacts.

### Thermal cycler parameters

Thermal cycling parameters were established for amplification of the SGM Plus™ Kit in the DNA Thermal Cycler 480 and GeneAmp™ PCR Systems 2400, 9600, and 9700. Thermal cycling times and temperatures met GeneAmp™ PCR Instrument specifications. Annealing and denaturation temperature windows were tested around each setpoint to verify that a  $\pm 2^\circ\text{C}$  window (DNA Thermal Cycler 480) or  $\pm 1.5^\circ\text{C}$  window (GeneAmp PCR System 2400, 9600, and 9700) yielded specific PCR product with the desired sensitivity of at least 1 ng of AmpF $\phi$ STR™ Control DNA 007.

SGM Plus™ Kit reactions were amplified for 27, 28, 29, 30, and 31 cycles on the GeneAmp PCR System 9600 using 1.5 ng of three DNA samples. As expected, PCR product increased with the number of cycles. Peak heights for all loci were above 150 RFU, even at 27 cycles; off-scale data was collected for several allele peaks at 31 cycles. See “DNA quantification” on page 19 for information on off-scale data.

While none of the cycle numbers tested produced nonspecific peaks, 28 cycles was found to give optimal sensitivity when the amplified products were examined on Applied Biosystems instruments. Additionally, the cycle number was set to avoid detection of low quantities of DNA (35 pg or less). At 28 cycles, 2.0 ng of AmpF $\phi$ STR™ Control DNA 007 amplifies reliably and specifically following the conditions outlined in this manual.

### AmpliTaq Gold™ DNA Polymerase activation

The thermal cycler program for the SGM Plus™ Kit includes an initial incubation to allow for AmpliTaq Gold™ DNA polymerase activation. Polymerase activation times and temperatures were tested in the DNA Thermal Cycler 480 and the GeneAmp PCR Systems 2400, 9600, and 9700.

Five activation times (5, 8, 11, 14 and 17 minutes) were tested. A plateau in fluorescent signal (RFU) of the SGM Plus™ Kit loci was reached at ~8 minutes and was maintained through the 17-minute timepoint. The 11-minute activation time was determined to be optimal based on this plateau (within a  $\pm >25\%$  window).

A 95°C activation temperature (for 11 minutes) was determined to be optimal for the reactions. These conditions produced maximum PCR product yield and a window in which reproducible performance characteristics were met.

## Accuracy, reproducibility, and precision

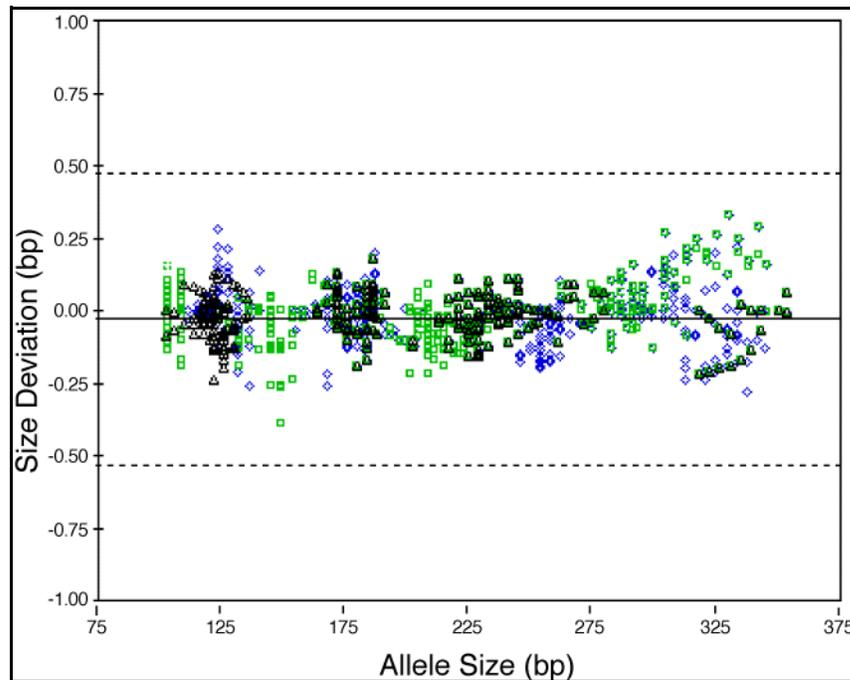
**DAB 8.1.2 Accuracy** *“Novel forensic DNA methodologies shall undergo developmental validation to ensure the accuracy, precision and reproducibility of the procedure.” (DAB, 1998).*

Laser-induced fluorescence detection systems of length polymorphism at short tandem repeat loci is not a novel methodology (Holt *et al.*, 2001 and Wallin *et al.*, 2001). However, accuracy and reproducibility of SGM Plus™ Kit profiles have been determined from various sample types.

## Accuracy

Figure 5 illustrates the size differences that are typically observed between sample alleles and AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder alleles on the Applied Biosystems 310 Genetic Analyzer with POP-4™ polymer. The x-axis in Figure 5 represents the nominal base pair sizes for a single injection of AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder, and the dashed lines parallel to the x-axis represent the  $\pm 0.5$ -bp windows. The y-axis is the deviation of each sample allele size from the corresponding allelic ladder allele size. The data include a total of 892 alleles from 30 population database samples and two injections of allelic ladder. In this representative example, all sample alleles are within 0.5 bp of a corresponding allelic ladder allele.

**Figure 5** Size deviation of 30 samples and two allelic ladders from one injection of allelic ladder on a single Applied Biosystems 310 instrument run.



The AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder contains the majority of alleles for the D3S1358, vWA, D16S539, D2S1338, amelogenin, D8S1179, D21S11, D18S51, D19S433, TH01, and FGA loci. However, alleles not found in the AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder do exist. These “off-ladder” alleles may contain full and/or partial repeat units. An “off-ladder” allele should flag itself by not falling inside the  $\pm 0.5$  bp window of any known allelic ladder allele.

**Note:** If a sample allele peak is found to be  $>0.5$  bp from the corresponding allelic ladder peak, then the sample must be rerun to verify the result.

## Reproducibility

The reproducible nature of purified DNA samples from various individuals ( $>10$ ), used routinely at Life Technologies, as well as validation samples processed during characterization of AmpF $\phi$ STR™ PCR Amplification Kits (for example, from body fluid mixture studies, environmental studies, matrix studies, nonprobative studies, CEPH family DNA sets) was without exception. All samples yielded the correct genotype.

## Precision and size windows

As indicated in the previous section, the recommended method for genotyping is to employ a  $\pm 0.5$ -bp “window” around the size obtained for each allele in the AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder. A  $\pm 0.5$ -bp window allows for the detection and correct assignment of potential off-ladder sample alleles whose true size is only one base different from an allelic ladder allele. Alleles of all possible sizes (within the range of 75–400 bp) should be readily identifiable. Any sample allele that sizes outside a window could be either of the following:

- An “off-ladder” allele, i.e., an allele of a size that is not represented in the AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder (go to <http://www.cstl.nist.gov/strbase/> for examples of known off-ladder alleles)
- An allele that does correspond to an allelic ladder allele, but whose size is just outside a window because of measurement error

The measurement error inherent in any sizing method can be defined by the degree of precision in sizing an allele multiple times. Precision is measured by calculating the standard deviation in the size values obtained for an allele that is run in several injections in one capillary run.

[Table 3 on page 70](#) indicates typical precision results obtained from 30 database samples and three AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder samples analyzed on the Applied Biosystems 310 Genetic Analyzer (47-cm capillary, POP-4™ polymer, GeneScan™ 500 ROX™ Size Standard). These results were obtained within a set of injections on a single capillary.

As indicated above, sample alleles may occasionally size outside of the  $\pm 0.5$ -bp window for a respective allelic ladder allele because of measurement error. The frequency of such an occurrence is lowest in detection systems with the smallest standard deviations in sizing. [Figure 5 on page 68](#) illustrates the tight clustering of allele sizes obtained on the Applied Biosystems 310 Genetic Analyzer, where the standard deviation in sizing is typically less than 0.15 bp. The instance of a sample allele sizing outside of the  $\pm 0.5$ -bp window because of measurement error is relatively rare when the standard deviation in sizing is approximately 0.15 bp or less (Smith, 1995).

For sample alleles that do not size within a  $\pm 0.5$ -bp window, the PCR product must be rerun to distinguish between a true off-ladder allele versus measurement error of a sample allele that corresponds with an allele in the allelic ladder.

It is important to note that while the precision within a set of capillary injections is very good, the determined allele sizes vary between platforms. Cross-platform sizing differences arise from a number of conditions, including type and concentration of polymer, run temperature, and electrophoresis conditions. Variations in sizing can also be found between runs on the same instrument and between runs on different instruments because of these conditions. We strongly recommend that the allele sizes obtained be compared to the sizes obtained for known alleles in the AmpF $\phi$ STR™ SGM Plus™ Allelic Ladder from the same run and then converted to genotypes (see [“Allelic ladder requirements” on page 25](#)). For more information on precision and genotyping, see Lazaruk *et al.*, 1998.

**Table 3** Example of precision results on the Applied Biosystems Genetic Analyzer

Allele	n	Mean	S.D.
<b>D3S1358</b>			
12	3	111.25	0.02
13	4	115.38	0.01
14	8	119.36	0.03
15	17	123.35	0.10
15	10	127.48	0.09
17	16	131.58	0.07
18	10	135.78	0.09
19	5	139.92	0.08
<b>vWA</b>			
11	3	154.59	0.04
12	3	158.80	0.01
13	3	163.00	0.02
14	11	167.22	0.11
15	7	171.18	0.05
16	13	175.15	0.07
17	12	179.18	0.06
18	16	183.15	0.06
19	9	187.09	0.07
20	3	190.94	0.04
21	3	194.88	0.05
22	3	198.73	0.06
23	3	202.55	0.06
24	3	206.83	0.04
<b>D16S539</b>			
5	3	229.29	0.06
8	4	241.27	0.01
9	8	245.25	0.08
10	6	249.33	0.07
11	19	253.27	0.07
12	22	257.31	0.06
13	12	261.36	0.07
14	3	265.51	0.04
15	3	269.57	0.03
<b>D2S1338</b>			
15	4	290.55	0.03
16	4	294.67	0.02

Allele	n	Mean	S.D.
17	11	298.75	0.06
18	8	303.12	0.07
19	11	307.53	0.07
20	10	311.86	0.10
21	4	316.20	0.05
22	7	320.39	0.12
23	9	324.68	0.15
24	11	328.72	0.09
25	10	332.93	0.11
26	4	336.92	0.12
27	3	340.89	0.06
28	3	344.57	0.08
<b>Amelogenin</b>			
X	29	103.48	0.06
Y	18	109.13	0.07
<b>D8S1179</b>			
8	3	123.69	0.02
9	6	127.73	0.03
10	7	131.80	0.08
11	6	135.92	0.40
12	9	140.20	0.06
13	19	144.72	0.10
14	14	149.23	0.10
15	14	153.70	0.07
16	3	158.08	0.02
17	3	162.27	0.06
18	3	166.38	0.04
19	3	170.48	0.04
<b>D21S11</b>			
24	3	185.21	0.03
24.2	3	187.21	0.04
25	3	189.17	0.03
26	3	193.08	0.05
27	3	197.01	0.06
28	11	200.82	0.07
28.2	3	202.78	0.06
29	12	204.77	0.05
29.2	3	206.71	0.04

<b>Allele</b>	<b>n</b>	<b>Mean</b>	<b>S.D.</b>
30	21	208.74	0.07
30.2	4	210.67	0.08
31	7	212.68	0.04
31.2	7	214.61	0.07
32	5	216.67	0.03
32.2	5	218.57	0.06
33	4	220.58	0.07
33.2	7	222.53	0.05
34	3	224.55	0.04
34.2	5	226.53	0.04
35	3	228.54	0.04
35.2	3	230.47	0.05
36	3	232.56	0.04
37	3	236.51	0.03
38	3	240.43	0.02
<b>D18S51</b>			
7	3	262.31	0.04
9	3	270.55	0.01
10	4	274.64	0.07
10.2	3	276.67	0.04
11	4	278.77	0.03
12	9	282.88	0.08
13	12	286.98	0.05
13.2	3	289.03	0.03
14	15	291.13	0.06
14.2	3	293.22	0.03
15	7	295.31	0.03
16	13	299.46	0.06
17	9	303.96	0.08
18	4	308.38	0.08
19	6	312.77	0.12
20	3	317.11	0.13
21	4	321.35	0.09
22	3	325.58	0.13
23	3	329.78	0.17
24	3	333.95	0.13
25	3	338.05	0.09
26	3	341.93	0.15

Allele	n	Mean	S.D.
27	3	345.59	0.10
<b>D19S433</b>			
9	3	102.01	0.05
10	3	105.93	0.04
11	3	109.82	0.07
12	6	113.68	0.06
12.2	3	115.73	0.04
13	16	117.62	0.04
13.2	4	119.65	0.06
14	22	121.57	0.08
14.2	6	123.60	0.07
15	14	125.54	0.06
15.2	5	127.62	0.06
16	7	129.54	0.08
16.2	3	131.59	0.06
17	3	133.60	0.05
17.2	3	135.62	0.03
<b>TH01</b>			
4	3	163.12	0.01
5	3	167.19	0.05
6	16	171.20	0.06
7	7	175.25	0.04
8	7	179.25	0.09
9	14	183.23	0.08
9.3	22	186.26	0.05
10	3	187.21	0.04
11	3	191.13	0.03
13.3	3	201.87	0.06
<b>FGA</b>			
17	3	212.48	0.02
18	5	216.51	0.07
19	7	220.52	0.06
20	10	224.54	0.06
21	16	228.52	0.07
22	14	232.60	0.07
23	12	236.62	0.03
24	8	240.64	0.07
25	8	244.74	0.06

Allele	n	Mean	S.D.
26	3	248.78	0.04
26.2	3	250.83	0.03
27	3	252.77	0.02
28	3	256.89	0.01
29	3	260.99	0.06
30	3	265.10	0.05
30.2	3	266.95	0.05
31.2	3	271.05	0.03
32.2	3	275.17	0.05
33.2	3	279.27	0.03
42.2	3	317.77	0.13
43.2	3	322.01	0.11
44.2	3	326.21	0.10
45.2	3	330.39	0.10
46.2	3	334.44	0.10
47.2	3	338.55	0.08
48.2	3	342.41	0.04
50.2	3	349.77	0.00
51.2	3	353.44	0.04

## Extra peaks in the electropherogram

### Overview

Peaks other than the target alleles may be detected on the electropherogram displays. Described below are several causes for the appearance of extra peaks, including the stutter product (found at the  $n-4$  position), incomplete 3' A nucleotide addition (found at the  $n-1$  position), and mixed DNA samples.

### Stutter products

The PCR amplification of tetranucleotide STR loci typically produces a minor product peak four bases shorter ( $n-4$ ) than the corresponding main allele peak. This is referred to as the stutter peak or product. Sequence analysis of stutter products at tetranucleotide STR loci has revealed that the stutter product is missing a single tetranucleotide core repeat unit relative to the main allele (Walsh *et al.*, 1996).

The proportion of the stutter product relative to the main allele (percent stutter) is measured by dividing the height of the stutter peak by the height of the main allele peak. Such measurements have been made for hundreds of samples at the loci used in the SGM Plus™ Kit.

Some of the general conclusions from these measurements and observations are as follows:

- For each SGM Plus™ Kit locus, the percent stutter generally increases with allele length, as shown in [Figure 6](#) through [Figure 8](#) on [page 76](#) through [page 76](#). Smaller alleles display a lower level of stutter relative to the longer alleles within each locus. This is reflected in [Figure 6](#) through [Figure 8](#), where minimal data points are plotted for some smaller alleles, because stutter was not detected for many of these samples.
- For the alleles within a particular locus, the percent stutter is generally greater for the longer allele in a heterozygous sample (this is related to the first point above).
- Each allele within a locus displays a percent stutter that is quite reproducible; the average of the standard deviation values measured for each allele at each locus is: 0.4% for TH01, 0.6% for D3S1358, vWA and FGA, 0.8% for D8S1179, D21S11 and D18S51, 0.9% for D16S539, and 1% for D19S433 and D2S1338. The expected range of percent stutter for any particular allele can be estimated as  $\pm 3$  standard deviations from the mean. For example, if the percent stutter for a particular allele averages 5% for multiple replicates, and if the average standard deviation at the allele is 0.5%, then the expected range in percent stutter for this allele is  $(5 \pm 1.5\%) = 3.5\text{--}6.5\%$ . This range also provides an estimate of the maximum expected stutter percent for each allele.
- The highest percent stutter observed for any TH01 allele was <4%, for any D3S1358, D8S1179, vWA or FGA allele <9%, for any D16S539 or D21S11 allele <10%, for any D2S1338 allele <12%, for any D18S51 allele <13%, and for any D19S433 allele <14%.
- An upper-limit stutter percent interpretational threshold can be estimated for each locus as 3 standard deviations above the highest percent stutter observed at the locus (see above two observations). Peaks at the stutter position that are above this threshold are not expected to be observed in single-source samples and therefore can be noted for closer examination. The upper-limit threshold values for each locus are as follows: 6% (TH01), 11% (D3S1358, vWA, FGA), 12% (D8S1179), 13% (D16S539, D21S11), 15% (D2S1338), 16% (D18S51), and 17% (D19S433). For evaluation of mixed samples, see [“DAB 8.1.2.2 Mixture Studies” on page 77](#).
- The percent stutter does not change significantly with the quantity of input DNA, for on-scale data with a minimum analyzed peak height of 150 relative fluorescence units (RFU). The measurement of percent stutter may be unnaturally high for main peaks that are off-scale. Loading or injecting less of the PCR product will yield accurate quantitation. See [“DNA quantification” on page 19](#) for information on off-scale data.

Figure 6 Stutter percentages for the D3S1358, vWA, D16S539, and D2S1338 loci

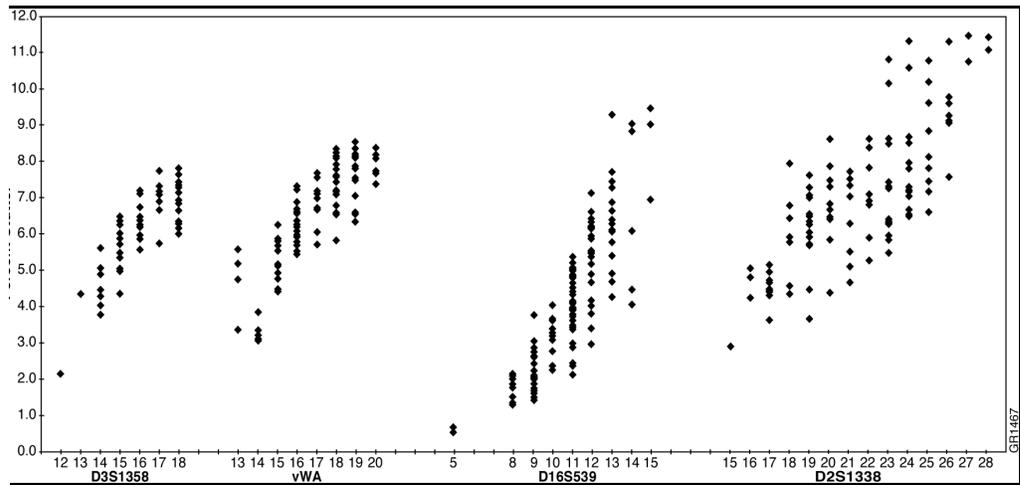


Figure 7 Stutter percentages for the D8S1179, D21S11, and D18S51 loci.

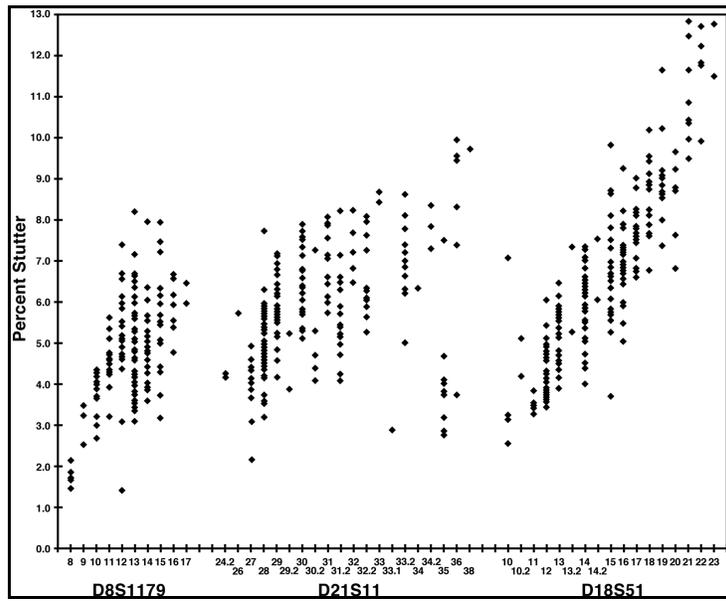
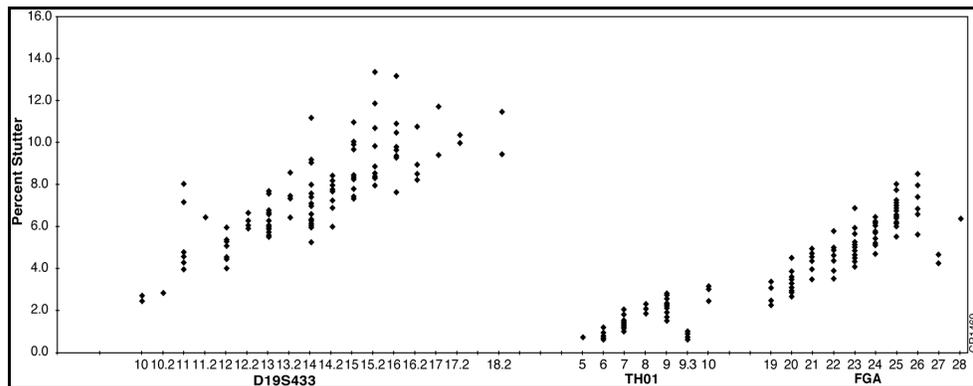


Figure 8 Stutter percentages for the D19S433, TH01, and FGA loci.



## Addition of 3' A nucleotide

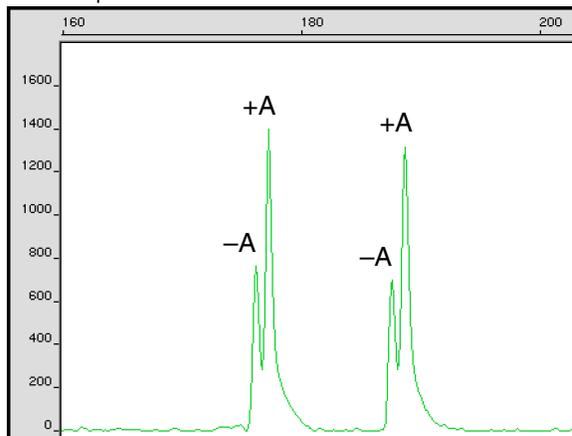
AmpliTaq Gold™ enzyme, like many other DNA polymerases, can catalyze the addition of a single nucleotide (predominately adenosine) to the 3' ends of double-stranded PCR products (Clark, 1988, Magnuson *et al.*, 1996). This non-template addition results in a PCR product that is one base pair longer than the actual target sequence, and the PCR product with the extra nucleotide is referred to as the “+A” form.

The efficiency of “A addition” is related to the particular sequence of the DNA at the 3' end of the PCR product. The SGM Plus™ Kit includes two main design features that promote maximum A addition:

- The primer sequences have been optimized to promote A addition.
- The last thermal cycling step is 60°C for 45 minutes.

This final extension step gives the AmpliTaq Gold™ DNA polymerase extra time to complete A addition to all double-stranded PCR product. STR systems that have not been optimized for maximum A addition may have “split peaks”, where each allele is represented by two peaks one base pair apart (Figure 9).

**Figure 9** Split peaks resulting from incomplete A nucleotide addition due to omission of the 45-minute extension step



Lack of full A nucleotide addition may be observed in SGM Plus™ Kit results when the amount of input DNA is greater than approximately 5.0 ng. The reason for this is that more time is needed for AmpliTaq Gold™ DNA Polymerase to add the A nucleotide to all molecules as more PCR product is generated. Amplification of too much input DNA will also result in off-scale data (see “DNA quantification” on page 19 for more information on off-scale data).

## DAB 8.1.2.2 Mixture Studies

“Species specificity, sensitivity, stability and mixture studies are conducted.” (DAB, 1998).

Evidence samples may contain DNA from more than one individual. The possibility of multiple contributors should be considered when interpreting the results. We recommend that individual laboratories assign a minimum peak height threshold based on validation experiments performed in each laboratory to avoid typing when stochastic effects are likely to interfere with accurate interpretation of mixtures.

In the discussion below, a peak is defined as any peak that is greater than 150 RFU. We recommend a minimum peak height threshold to avoid typing less than 35 pg of input DNA (see “Effect of DNA quantity on results” on page 85).

## Detection of mixed samples

Each of the following can aid in determining whether a sample is a mixture:

- The presence of more than two alleles at a locus
- The presence of a peak at a stutter position that is significantly greater in percentage than what is typically observed in a single-source sample (see “Stutter products” on page 74 and Figure 6 through Figure 8).
- Significantly imbalanced alleles for a heterozygous genotype  
The peak height ratio is defined as the height of the lower peak (in RFU) divided by the height of the higher peak (in RFU), expressed as a percentage. Mean peak height ratios and standard deviations observed for alleles in the SGM Plus™ Kit loci in unmixed population database samples are as follows:

Allele	Mean Peak Height Ratio	Number of Observations (n)
D3S1358	93 ± 4%	68
vWA	93 ± 5%	74
D16S539	92 ± 6%	70
D2S1338	91 ± 6%	87
Amelogenin	90 ± 6%	46
D8S1179	92 ± 6%	93
D21S11	91 ± 7%	95
D18S51	91 ± 6%	100
D19S433	92 ± 6%	97
TH01	92 ± 6%	70
FGA	93 ± 5%	80

- For all 11 loci, the mean peak height ratios indicate that the two alleles of a heterozygous individual are generally very well balanced. Ratios <70% are rare in normal, unmixed samples.

If the peak height ratio is <70% for one locus, and there are no other indications that the sample is a mixture, the sample may be reamplified and reanalyzed to determine if the imbalance is reproducible. Reproducible imbalance at only one locus may indicate a mixture of significantly overlapping genotypes. Other possible causes of imbalance at a locus are degraded DNA, presence of inhibitors, extremely low amounts of input DNA, or the presence of an allele containing a rare sequence that does not amplify as efficiently as the other allele. Amplification and analysis of additional loci may assist in the interpretation of the sample.

## Resolution of genotypes in mixed samples

A sample containing DNA from two sources can be comprised (at a single locus) of any of the seven genotype combinations listed below.

- Heterozygote + heterozygote, no overlapping alleles (four peaks)
- Heterozygote + heterozygote, one overlapping allele (three peaks)
- Heterozygote + heterozygote, two overlapping alleles (two peaks)
- Heterozygote + homozygote, no overlapping alleles (three peaks)
- Heterozygote + homozygote, overlapping allele (two peaks)
- Homozygote + homozygote, no overlapping alleles (two peaks)
- Homozygote + homozygote, overlapping allele (one peak)

Specific genotype combinations and input DNA ratios of the samples contained in a mixture determine whether it is possible to resolve the genotypes of the major and minor component(s) at a single locus.

The ability to obtain and compare quantitative values for the different allele peak heights on Life Technologies instruments provides additional valuable data to aid in resolving mixed genotypes. This quantitative value is much less subjective than comparing relative intensities of bands on a stained gel.

Ultimately, the likelihood that any sample is a mixture must be determined by the analyst in the context of each particular case, including the information provided from known reference sample(s).

## Limit of detection of the minor component

Mixtures of two DNA samples were examined at various ratios (1:1 to 1:20). The total amount of genomic input DNA mixed at each ratio was 1 ng.

The samples were amplified in a GeneAmp™ PCR System 9600 and were electrophoresed and detected using a 377 DNA Sequencer.

The results of the mixed DNA samples, shown separately in [Figure 10 on page 80](#), are displayed in [Figure 11 on page 81](#) where sample A was the minor component and sample B was the major component.

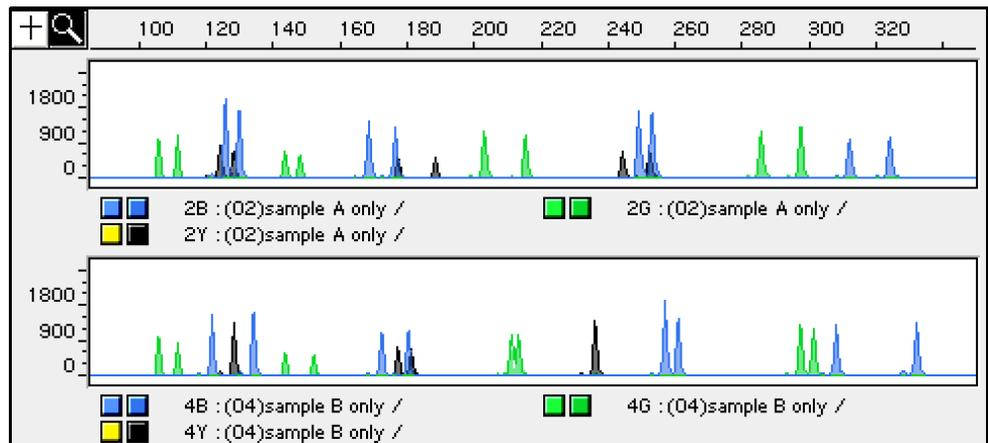
The genotypes of the samples in [Figure 10](#) are the following:

Allele	Genotype	
	Sample A	Sample B
Amelogenin	X, Y	X, Y
D19S433	14,15	15,15
D3S1358	15, 16	14,17
D8S1179	12, 13	12,14
vWA	14, 16	15,17
TH01	7, 9.3	7, 8
D21S11	28, 31	30, 30.2
FGA	24, 26	22, 22
D16S539	9, 10	11, 12
D18S51	12, 15	15, 16
D2S1338	20, 23	19, 25

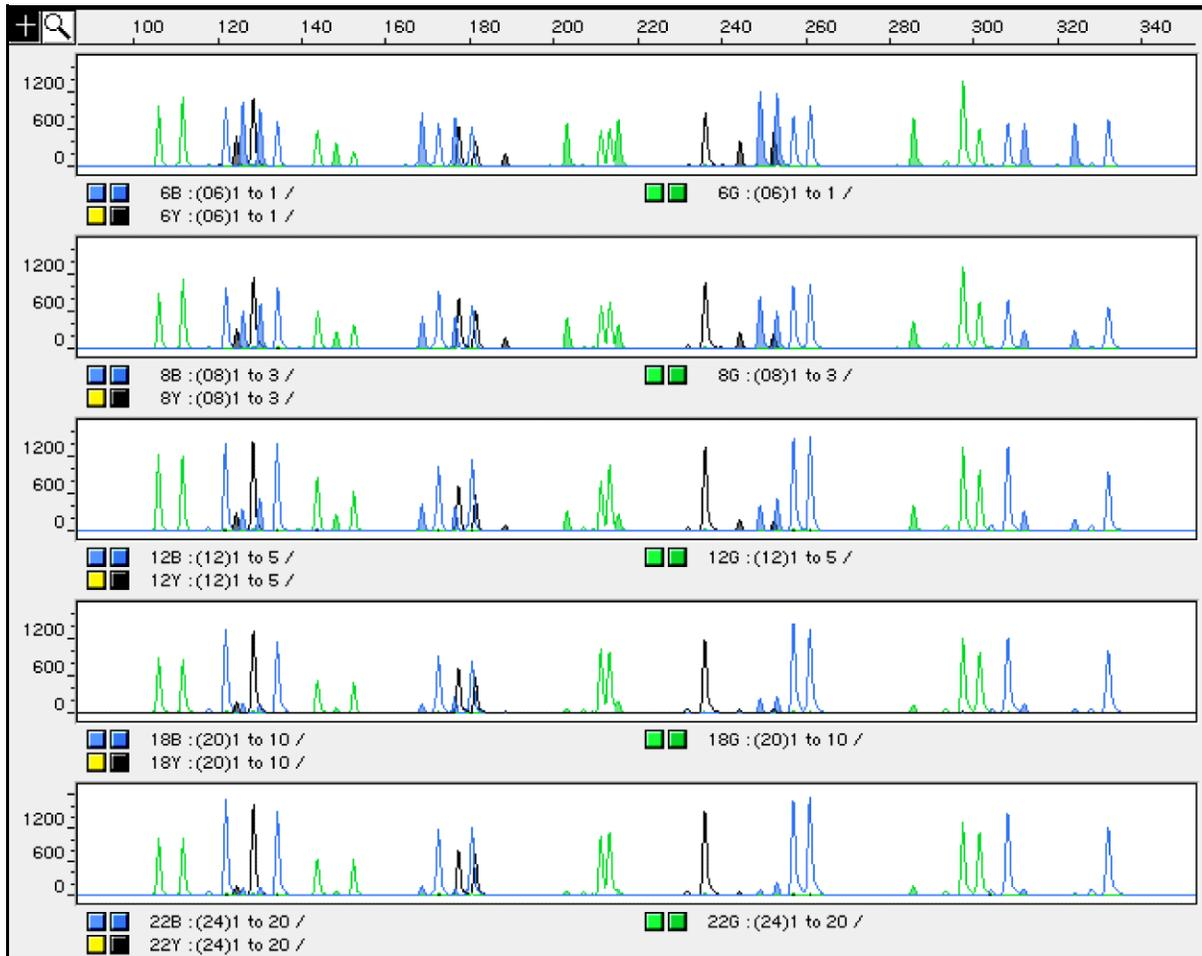
For these 1 ng total DNA mixture studies, the limit of detection is when the minor component is present at approximately 1/20 of the concentration of the major component. The limit of detection for the minor component is influenced by the combination of genotypes in the mixture.

The following figures show the reference samples and the two DNA samples used for this study.

**Figure 10** Reference samples for mixture study shown in next figure



**Figure 11** Results of the two DNA samples from previous figure mixed together at defined ratios and amplified with the SGM Plus™ Kit. The A:B ratios shown are 1:1, 1:3, 1:5, 1:10, and 1:20 (top to bottom). Alleles attributable only to the minor component are highlighted.



Simulation of forensic casework scenarios was achieved by combining various body fluids (blood: blood, semen: blood, and semen: saliva) from two donors in defined ratios, by volume, from 1:1–1:50.

Mixed stains were prepared and DNA was extracted at the Santa Clara County Crime Laboratory, San Jose, CA, following the phenol/chloroform procedure (differential lysis was performed on stains containing semen). At Applied Biosystems approximately 3 ng of DNA was amplified from each dried fluid mixture in a GeneAmp PCR System 9600 and detected using either the Applied Biosystems 310 Genetic Analyzer or the Applied Biosystems 377 DNA Sequencer.

The limit of detection of the minor genotype component of each body fluid mixture ratio was determined. The limit of detection is defined here as the ratio below which a mixture is not recognized but rather appears to contain DNA from a single source (the major contributor of the mixture). The limit of detection occurred in blood: blood mixtures when the minor component was present at one tenth the volume of the major genotype.

In epithelial cell fractions of differential extractions of semen: blood or semen: saliva stains, blood was detectable at one fiftieth and saliva at one tenth the volume of semen. Sperm DNA carryover into the epithelial cell fraction of these stains was detectable at one-tenth the volume of blood or saliva. In sperm fractions, the male genotype was detectable from every semen: blood or semen: saliva mixture with no trace of the female DNA.

Again, the limit of detection for the minor component is influenced by the specific combination of genotypes present in mixtures.

Note that the limit of detection for the minor component is influenced by the specific combination of genotypes present in mixtures.

### Analysis of sexual assault DNA mixture evidence

SGM Plus™ Kit reactions with DNA extracted from adjudicated sexual assault evidence were examined to assess the performance of the SGM Plus™ Kit on typical casework samples comprised of mixed body fluids. These samples were extracted, amplified, and analyzed in collaboration with the Santa Clara County Crime Laboratory.

DNA extracts from three adjudicated sexual assault cases were prepared by DNA analysts at the Santa Clara County Crime Laboratory. Sexual assault evidence materials were processed using the differential lysis and organic extraction procedure, while victim/suspect reference blood samples were processed using the Chelex™ extraction procedure. Following amplification with the SGM Plus™ Kit reagents, the PCR products were analyzed using the Applied Biosystems 377 DNA Sequencer.

- Case 1 and Case 2 contained a victim reference blood sample, a suspect reference blood sample, and a victim vaginal swab. The SGM Plus™ Kit genotype of the epithelial cell fraction was the same as that of the victim reference and did not contain alleles foreign to the victim. The SGM Plus™ Kit genotype of the sperm cell fraction did not contain detectable epithelial cell fraction DNA and included the suspect as a possible semen donor.
- Case 3 contained a victim reference blood sample and a victim vaginal swab. The SGM Plus™ Kit genotype of the epithelial cell fraction was the same as that of the victim reference and did not contain alleles foreign to the victim. In accordance with the victim's account, SGM Plus™ Kit genotypes of the sperm fraction revealed DNA from multiple semen donors. No suspect(s) were developed in this case.

## Characterization of loci

### DAB 8.1.2.1 Documentation

*"Documentation exists and is available which defines and characterizes the locus."* (DAB, 1998).

### Nature of the polymorphisms

The primers for the amelogenin locus flank a six-base pair deletion within intron 1 of the X homologue. Amplification results in 107-bp and 113-bp products from the X and Y chromosomes, respectively. (Sizes are the actual base pair size according to sequencing results, including 3' A nucleotide addition.) The remaining SGM Plus™ Kit loci are all tetranucleotide short tandem repeat (STR) loci. The length differences among alleles of a particular locus result from differences in the number of 4-bp repeat units (see [Table 1 on page 12](#)).

Alleles in the AmpFtSTR™ SGM Plus™ Allelic Ladder, alleles containing partial repeat units, population database, and nonhuman primate DNA samples have been subjected to DNA sequencing at Life Technologies. In addition, other groups in the forensic community have sequenced alleles at some of these loci (Nakahori *et al.*, 1991; Puers *et al.*, 1993; Möller *et al.*, 1994; Barber *et al.*, 1995; Möller and Brinkmann, 1995; Barber *et al.*, 1996; Barber and Parkin, 1996; Brinkmann *et al.*, 1998; Momhinweg *et al.*, 1998; Watson *et al.*, 1998). Among the various sources of sequence data on the SGM Plus™ Kit loci, there is consensus on the repeat patterns and structure of the STRs (see [Table 1 on page 12](#)).

D19S433 alleles have been sequenced at Life Technologies, including those in the AmpFtSTR™ SGM Plus™ Allelic Ladder. Sequenced D19S433 alleles that contain a 2-bp repeat unit result from a deletion of the last two nucleotides (AG) in the second tetranucleotide repeat unit (AAAG) of the repeat motif (see [Table 1 on page 12](#)).

## Inheritance

The SGM Plus™ Kit loci have been validated by family studies in order to demonstrate their mode(s) of inheritance.

The Centre d'Etude du Polymorphisme Humain (CEPH) has collected DNA from 39 families of Utah Mormon, French Venezuelan, and Amish descent. These DNA sets have been extensively studied all over the world and are routinely used to characterize the mode of inheritance of various DNA loci. Each family set contains three generations, generally including four grandparents, two parents, and several offspring. Consequently, the CEPH family DNA sets are ideal for studying inheritance patterns (Begovich *et al.*, 1992).

Four CEPH family DNA sets were examined. DNA input of 1.5 ng from each sample was amplified using the SGM Plus™ Kit, followed by analysis using an Applied Biosystems 377 DNA Sequencer. The families examined included #1331 (11 offspring), #13291 (9 offspring), #13292 (9 offspring), and #13294 (8 offspring), representing 37 meiotic divisions. The results confirmed that the loci are inherited according to Mendelian rules, as has been reported in the literature (Nakahori *et al.*, 1991; Edwards *et al.*, 1992; Kimpton *et al.*, 1992; Mills *et al.*, 1992; Sharma and Litt, 1992; Li *et al.*, 1993; Straub *et al.*, 1993).

## Mapping

The SGM Plus™ Kit loci D3S1358, vWA, D16S539, D2S1338, amelogenin, D8S1179, D21S11, D18S51, D19S433, TH01, and FGA have been mapped and the chromosomal locations have been published (Nakahori *et al.*, 1991; Edwards *et al.*, 1992; Kimpton *et al.*, 1992; Mills *et al.*, 1992; Sharma and Litt, 1992; Li *et al.*, 1993; Straub *et al.*, 1993; Barber and Parkin, 1996). They are listed in [Table 1 on page 12](#).

## Population studies

Population distribution data of the 11 SGM Plus™ Kit loci have been established in different racial and/or ethnic groups. These loci were amplified and typed for 200 U.S. Caucasian and 195 African-American individuals. For more information regarding analysis of these samples, see ["Population data" on page 92](#)

## Species specificity

### DAB 8.1.2.2 Species Specificity

*"Species specificity, sensitivity, stability and mixture studies are conducted."* (DAB, 1998).

**Nonhuman studies**

Nonhuman DNA may be present in forensic casework samples. The SGM Plus™ Kit provides the required degree of specificity such that it is specific to primates for the species tested (with the exception of the amelogenin locus).

The following experiments were conducted to investigate interpretation of SGM Plus™ Kit results from nonhuman DNA sources.

The extracted DNA samples were amplified in SGM Plus™ Kit reactions and analyzed using the Applied Biosystems 377 DNA Sequencer.

- **Primates** – Gorilla, chimpanzee, and macaque (2.5 ng each).
- **Non-primates** – Hamster, rat, rabbit, cat, dog, pig, chicken, cow, and fish (50 ng each).
- **Bacteria and yeast** – *Legionella*, *Escherichia*, *Listeria*, *Neisseria*, *Vibrio*, *Citrobacter*, *Salmonella*, *Candida*, *Saccharomyces* (equivalent to ~50 ng human DNA), and *Rhodotorula*.

The primate DNA samples all amplified, producing fragments within the 75–350 base pair region (Wallin *et al.*, 1998). The primate samples were subsequently sequenced by Life Technologies scientists. The data revealed significant sequence homology between the primate and human DNA for the SGM Plus™ Kit loci (Meyer *et al.*, 1995; Andersen *et al.*, 1996).

The bacteria, yeast, hamster, rat, rabbit, cat, chicken, and fish samples did not yield detectable product. The dog, pig, and cow samples produced a 103-bp fragment. This 102-bp fragment was also amplified using the amelogenin primers alone. This confirms amplification of the product obtained by Buel *et al.* (1995). The 103-bp fragment is 4 bp shorter than the primate 107-bp X-specific product (including +A addition).

## Sensitivity

### DAB 8.1.2.2 Sensitivity

*“Species specificity, sensitivity, stability and mixture studies are conducted.”* (DAB, 1998).

### Effect of DNA quantity on results

The amount of input DNA added to the PCR reaction should be 1.0–2.5 ng. The DNA sample should be quantitated prior to amplification using a system such as the Quantifiler™ Human DNA Quantitation Kit (Part no. 4343895). [Figure 12 on page 86](#) shows the effect of different amounts of AmpFtSTR™ Control DNA 007.

The final DNA concentration should be in the range of 0.05–0.125 ng/μL so that 1.0–2.5 ng of DNA will be added to the PCR reaction in a volume of 20 μL. If the sample contains degraded DNA, amplification of additional DNA may be beneficial.

If too much DNA is added to the PCR reaction, then the increased amount of PCR product that is generated can result in the following:

- Fluorescence intensity that exceeds the linear dynamic range for detection by the instrument (“off-scale” data)

Off-scale data is a problem for two reasons:

- Quantitation (peak height and area) for off-scale peaks is not accurate. For example, an allele peak that is off-scale can cause the corresponding stutter peak to appear higher in relative intensity, thus increasing the calculated percent stutter.
- Multicomponent analysis of off-scale data is not accurate, which results in poor spectral separation (“pull-up”).

Identification of off-scale peaks and multicomponent analysis are discussed in [“DNA quantification” on page 19](#) and [“About multicomponent analysis” on page 16](#).

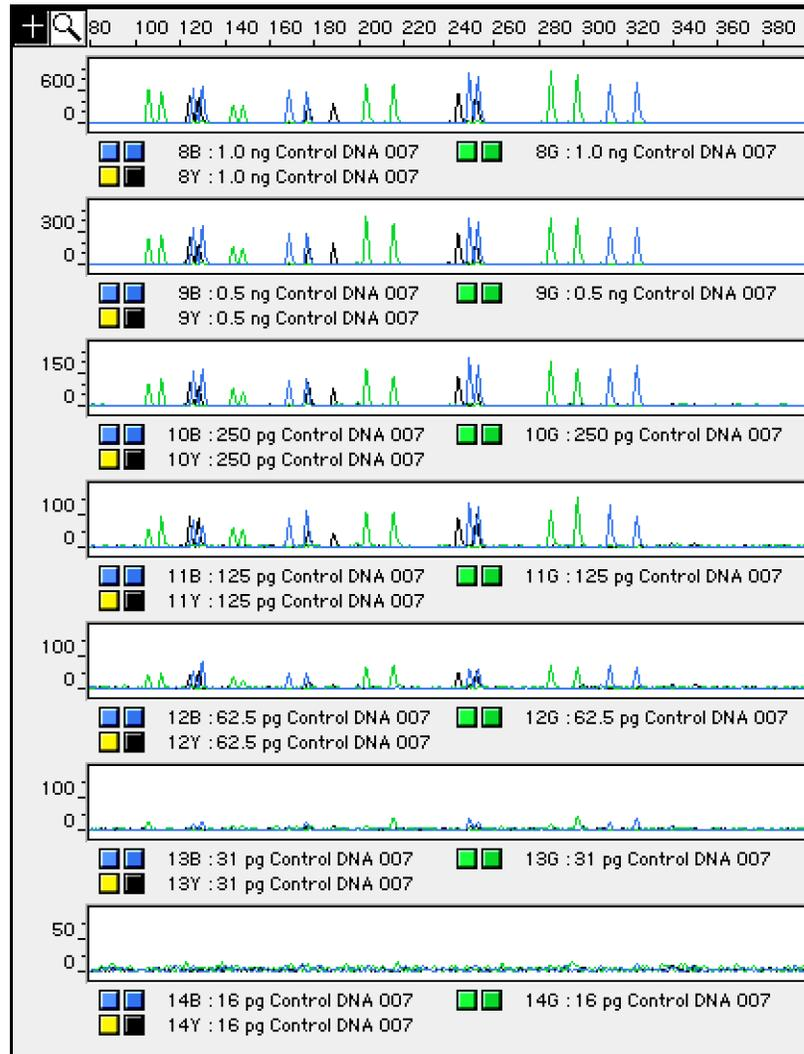
- Incomplete A nucleotide addition

To avoid these issues, the sample can be re-amplified using less DNA.

When the total number of allele copies added to the PCR is extremely low, unbalanced amplification of the two alleles of a heterozygous individual may occur (Wallin *et al.*, 1998; Walsh *et al.*, 1992). This is due to stochastic fluctuation in the ratio of the two different alleles (Sensabaugh *et al.*, 1991). The PCR cycle number and amplification conditions have been specified to produce low peak heights for a sample containing produce peak heights of <150 RFU for a sample containing 35 pg human genomic DNA (corresponding to ten total allele copies). Peak heights <150 RFU should be interpreted with caution.

Individual laboratories may find it useful to determine an appropriate minimum peak height interpretational threshold based on their own results using low amounts of input DNA.

**Figure 12** Effect of amplifying various amounts of AmpF $\phi$ STR™ Control DNA 007 ranging from 16 pg to 1 ng. Note that the y-axis scale differs in many of these panels.



## Stability

### DAB 8.1.2.2 Stability

*"Species specificity, sensitivity, stability and mixture studies are conducted."* (DAB, 1998).

### Lack of amplification of some loci

As with any multi-locus system, the possibility exists that not every locus will amplify. This is most often observed when the DNA substrate has been severely degraded or when the DNA sample contains PCR inhibitors. Because each locus is an independent marker, results generally can still be obtained from the loci that do amplify.

## Differential and preferential amplification

Differential amplification can be defined as the difference in the degree of amplification of each locus within a co-amplified system, such that one or more loci may amplify to a greater extent compared to the other loci. Preferential amplification is used in this guide to describe differences in the amplification efficiency of two alleles at a single locus and is observed when the peak height ratio between the two alleles is <70% (see [“DAB 8.1.2.2 Mixture Studies” on page 77](#)).

Preferential amplification of alleles in systems that distinguish alleles based on length polymorphisms is most likely to be observed when the alleles differ significantly in base pair size. Because most SGM Plus™ Kit loci have small size ranges, the potential for preferential amplification of alleles is low.

DNA samples containing either D2S1338 alleles separated by 28–40 bp or containing FGA alleles separated by 90–100 bp were analyzed in several of our studies to determine the potential for preferential amplification.

In assessing potential for differential and preferential amplification, the following four variables were examined:

- Low template copy number
- Effect of PCR inhibitor in a DNA sample
- Degraded DNA
- Amplification denaturation and annealing temperatures

### Low template copy number

To determine if the amount of input DNA may result in either differential or preferential amplification, varying quantities of three DNA samples were amplified in duplicate reactions. One nanogram, 0.5 ng, 0.25 ng, 0.125 ng, 0.06 ng, 0.03 ng, and 0.015 ng of AmpF $\phi$ STR Control DNA 007 and samples containing either widespread D2S1338 or FGA alleles were amplified and then analyzed using the Applied Biosystems 377 DNA Sequencer.

No loci in any of the samples tested differentially amplified at any of the eleven loci. Preferential amplification was not observed between the widespread D2S1338 alleles.

For the sample containing widespread FGA alleles, the duplicate dilution series are denoted here as A and B. In dilution series A, for the sample containing 0.06 ng template DNA, the shorter of the two widespread FGA alleles (270 bp) was detected and the longer allele (334 bp) was not. However, the shorter allele peak height (26 RFU) was significantly below the recommended minimum threshold of 150 RFU (see [“Effect of DNA quantity on results” on page 85](#)). In dilution series A, preferential amplification was not seen at any of the higher template DNA concentrations.

Moreover, in dilution series B, for the sample containing widespread FGA alleles, such preferential amplification did not occur at 0.06 ng. But, at 0.125 ng of input DNA, a peak height ratio of 55% was detected between the widespread FGA alleles. Both FGA alleles in this sample (93 and 51 RFU) were below the recommended minimum threshold of 150 RFU (see [“Effect of DNA quantity on results” on page 85](#)).

For dilution series A and B, at 0.25 ng and above, both FGA alleles amplified in equal proportions; at 0.03 ng neither allele was detectable. Taken together, these data underscore the importance of interpreting single peaks of low fluorescence signal with caution and on a case-by-case basis.

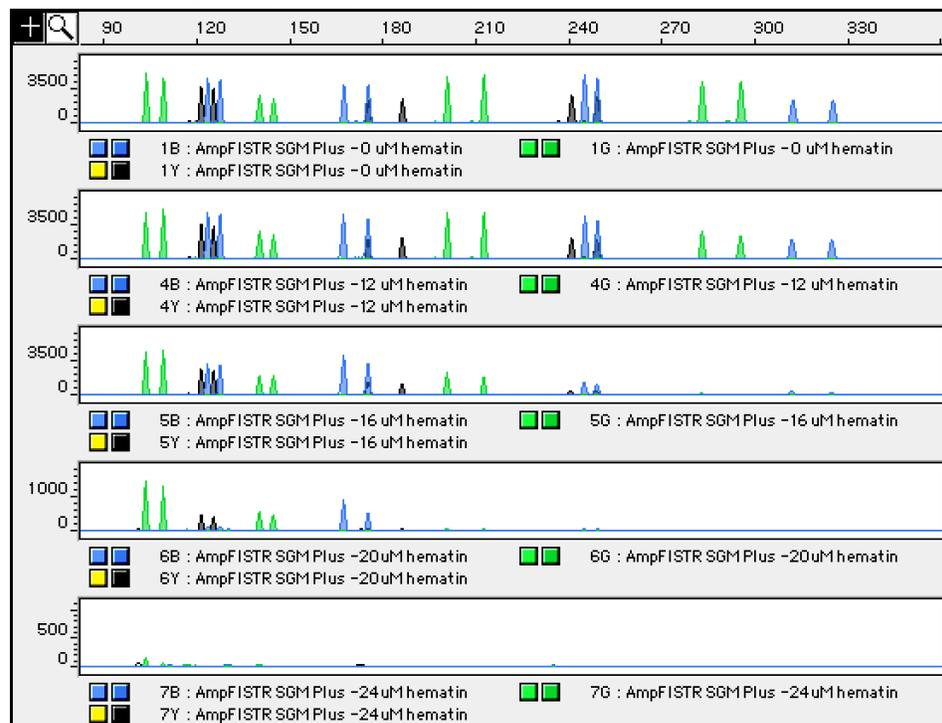
## Effect of inhibitors

Heme compounds have been identified as PCR inhibitors in DNA samples extracted from bloodstains (Akane *et al.*, 1994., DeFranchis *et al.*, 1988). It is believed that the inhibitor is co-extracted and co-purified with the DNA and subsequently interferes with PCR by inhibiting polymerase activity.

Bovine serum albumin (BSA) can prevent or minimize the inhibition of PCR, most likely by binding to the inhibitor (Comey *et al.*, 1994). Since the presence of BSA can improve the amplification of DNA from blood-containing samples, BSA has been included in the AmpF $\phi$ STR™ PCR Reaction Mix at a concentration of 8  $\mu$ g per 50- $\mu$ L amplification. BSA has also been identified as an aid in overcoming inhibition from samples containing dyes, such as in denim.

To examine the effects of hematin on the SGM Plus™ Kit amplification results, DNA samples were amplified using the SGM Plus™ Kit reagents (including the BSA-containing PCR reaction mix) in the presence of varying concentrations of purified hematin. The concentrations of hematin used were 0  $\mu$ M, 12  $\mu$ M, 16  $\mu$ M, 20  $\mu$ M, and 24  $\mu$ M. When the amount of hematin was increased to a concentration that started to inhibit the PCR, D2S1338 and D18S51 were the first loci to drop out in each experiment, followed by FGA (Figure 13).

**Figure 13** DNA amplified with the SGM Plus™ Kit in the presence of varying concentrations of hematin: 0, 12  $\mu$ M, 16  $\mu$ M, 20  $\mu$ M, and 24  $\mu$ M

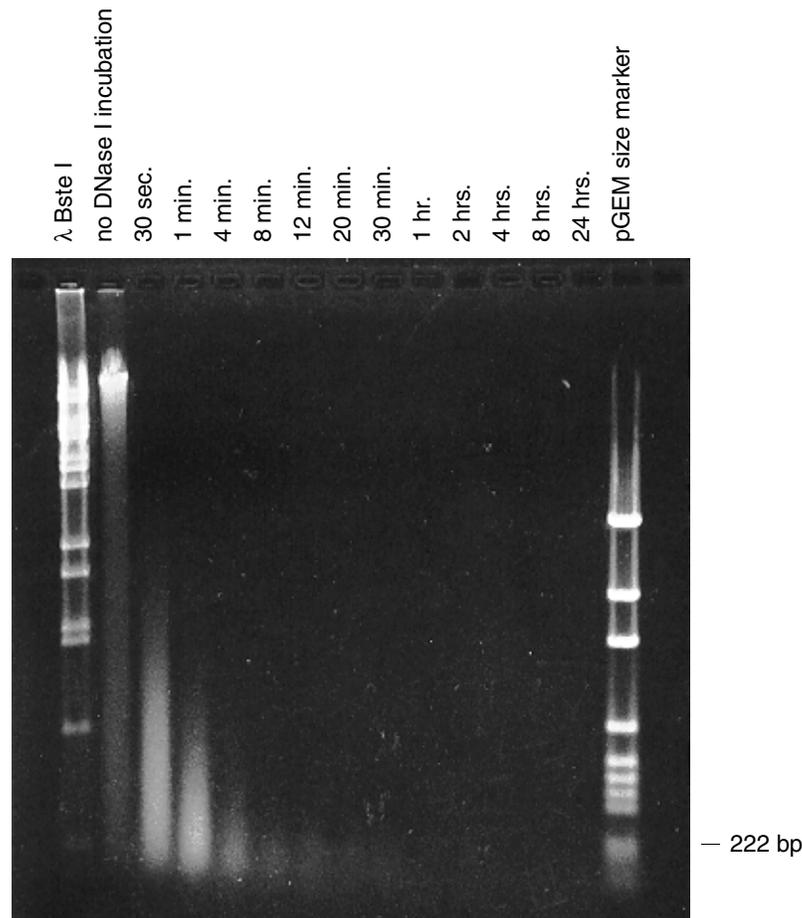


## Degraded DNA

As the average size of degraded DNA approaches the size of the target sequence, the amount of PCR product generated is reduced. This is due to the reduced number of intact templates in the size range necessary for amplification.

Degraded DNA was prepared to examine the potential for differential amplification of loci. High molecular weight DNA was incubated with the enzyme DNase I for varying amounts of time. The DNA was examined by agarose gel analysis to determine the average size of the DNA fragments at each timepoint. Gel analysis results of the degraded DNA are shown in [Figure 14](#).

**Figure 14** Agarose gel of degraded genomic DNA



Four (4) ng of degraded DNA (or 2 ng undegraded DNA) was amplified using the SGM Plus™ Kit (all 11 primer pairs together) and also in reactions containing each locus-specific primer pair individually.

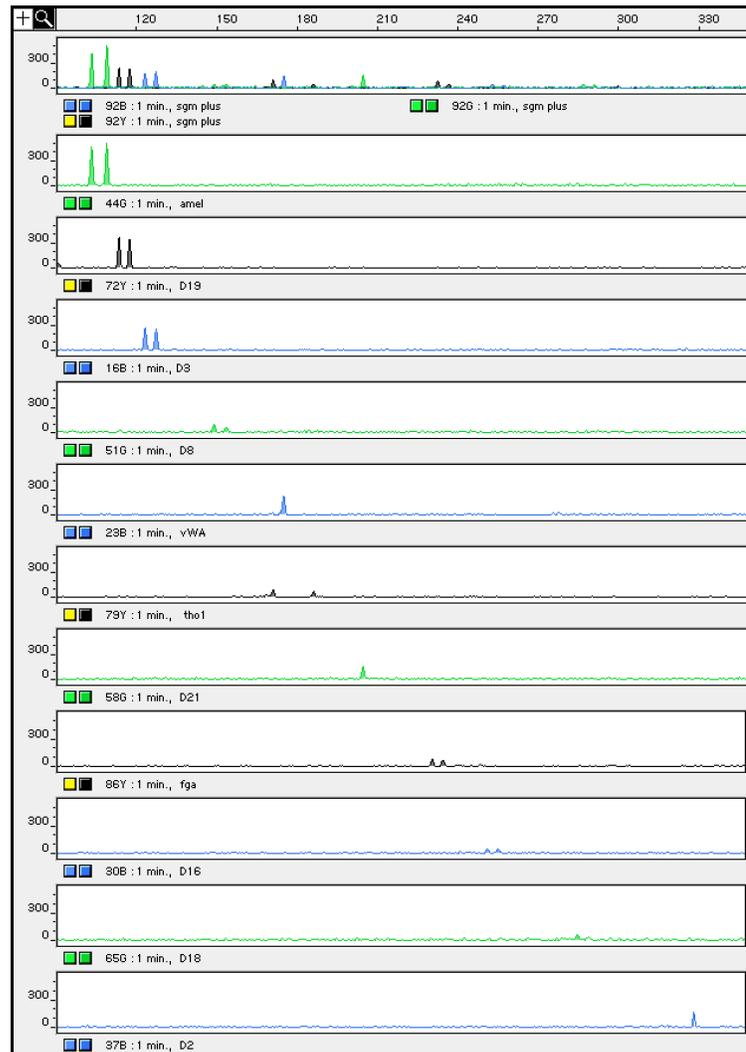
The electropherograms in [Figure 15](#) show the SGM Plus™ Kit amplification results of a DNA sample with no DNase I treatment (2 ng amplified) and those of the 30-second, 1-, 4-, and 8-minute incubations (approximately 4 ng amplified).

Figure 15 Amplifications of DNA incubated for various times with DNase I



The loci failed to amplify in the order of decreasing size as the extent of degradation progressed: D2S1338 was the first locus to drop out, followed by D18S51 and D16S539 and so forth. A similar result at each timepoint was obtained whether the DNA samples were amplified for each locus alone or co-amplified with the SGM Plus™ Kit (Figure 16).

**Figure 16** Multiplex and single-locus amplifications of the DNA sample incubated for 1 minute with DNase I



### Denaturation and annealing temperatures

The effects of denaturation and annealing temperatures on the amplification of SGM Plus™ Kit loci were examined using 1–2 ng of three DNA samples: AmpFtSTR Control DNA 007 and samples containing either widespread D2S1338 or FGA alleles.

The denaturation temperatures tested were 92, 94, and 96°C (DNA Thermal Cycler 480) and 92.5, 94, and 95.5 °C (GeneAmp Detection Systems 2400, 9600, and 9700), all for 1-minute hold times. The annealing temperatures tested were 57, 59, 61, and 63°C, also for 1-minute hold times, in the DNA Thermal Cycler 480 and the GeneAmp PCR Systems 2400, 9600 and 9700. The PCR products were analyzed using the Applied Biosystems 310 Genetic Analyzer.

Neither preferential nor differential amplification was observed in the denaturation temperature experiments. Of the tested annealing temperatures, 57, 59, and 61°C did not induce any differential amplification. At 63°C, the yield of the majority of loci was significantly reduced. This should pose no problem with routine thermal cyclers calibration and when following the recommended amplification protocol. Preferential amplification was not observed at any of the tested annealing temperatures.

## Matrix studies

Analysts at the Santa Clara Crime Laboratory prepared a panel of blood and semen specimens deposited on a variety of commonly encountered substrates. Blood samples from two donors were deposited individually on wool, cotton, nylon, metal, glass, leather (one donor), and blue denim. Semen samples from two donors were deposited individually on wool, cotton, nylon, leather, blue denim, acetate, vinyl upholstery, facial tissue, and a condom with spermicide (5% nonoxynol-9).

Specimens were stored at room temperature and at specified time points a sampling of each stain was removed for extraction. The blood and semen stains were extracted using the organic extraction procedure. Additionally, a portion of each blood specimen was extracted using the Chelex™ method followed by Centricon™-100 ultrafiltration. Extracted samples were then stored at –15 to –25°C for 4–5 years.

The 1–week and 1–year time points were analyzed by Life Technologies scientists. The samples were amplified using SGM Plus™ Kit reagents and analyzed using the Applied Biosystems 310 Genetic Analyzer. A complete 11-locus genotype was obtained for 47 of the 49 blood and semen samples exposed to the tested matrices for 1 week or 1 year. Some alleles in two of the 1–year samples were of low signal intensity and were either too weak to be genotyped or were interpreted with caution. No non-specific artifacts were generated and all samples yielded expected genotypes.

## Population data

### DAB 8.1.2.3 Population Data

*“Population distribution data are documented and available.” (DAB, 1998).*

### DAB 8.1.2.3.1 Population Distribution Data

*“The population distribution data would include the allele and genotype distributions for the locus or loci obtained from relevant populations. Where appropriate, databases should be tested for independence expectations.” (DAB, 1998).*

### Overview

To interpret the significance of a match between genetically typed samples, it is necessary to know the population distribution of alleles at each locus in question. If the genotype of the relevant evidence sample is different from the genotype of the suspect's reference sample, then the suspect is “excluded” as the donor of the biological evidence tested. An exclusion is independent of the frequency of the two genotypes in the population.

If the suspect and evidence samples have the same genotype, then the suspect is “included” as a possible source of the evidence sample. The probability that another, unrelated, individual would also match the evidence sample is estimated by the frequency of that genotype in the relevant population(s).

**Population samples used in these studies**

The SGM Plus™ Kit was used to generate the population data provided in this section. Samples were collected from individuals throughout the United States with no geographical preference.

Population	Number of samples	Samples provided by
African-American	195	Laboratory Corporation of America
U.S. Caucasian	200	

**Allele frequencies**

Table 4 shows the SGM Plus™ Kit allele frequencies in two populations, listed as percentages.

**Table 4** SGM Plus™ Kit allele frequencies

Allele	African-American (n = 195)	U.S. Caucasian (n = 200)
<b>D3S1358</b>		
9	0.26 <sup>†</sup>	‡
11	0.26 <sup>†</sup>	0.25 <sup>†</sup>
12	0.51 <sup>†</sup>	0.25 <sup>†</sup>
13	‡	0.50 <sup>†</sup>
14	11.80	11.25
15	27.95	28.25
15.2	0.26 <sup>†</sup>	‡
16	32.31	22.25
17	21.80	22.25
18	4.62	14.50
19	0.26 <sup>†</sup>	0.50 <sup>†</sup>
<b>vWA</b>		
11	0.26 <sup>†</sup>	‡
12	‡	‡
13	1.54	‡
14	7.70	8.50
15	22.05	8.25
16	26.92	19.75
17	16.92	25.00
18	13.85	25.75
19	8.46	11.00
20	2.05	1.50
21	‡	0.25 <sup>†</sup>
22	0.26 <sup>†</sup>	‡
23	‡	‡
24	‡	‡

Allele	African-American (n = 195)	U.S. Caucasian (n = 200)
<b>D16S539</b>		
5	0.26 <sup>†</sup>	‡
8	3.85	1.5
9	18.21	10.75
10	11.28	4.25
11	30.00	29.75
12	18.46	34.25
13	15.64	17.5
14	2.05	1.75
15	0.26 <sup>†</sup>	0.25 <sup>†</sup>
<b>D2S1338</b>		
15	‡	0.25 <sup>†</sup>
16	5.64	4.25
17	11.03	17.75
18	2.82	8.75
19	16.41	12.75
20	8.72	13.50
21	13.59	3.25
22	12.31	3.75
23	8.72	10.75
24	8.46	10.00
25	8.72	12.00
26	2.82	2.75
27	0.51 <sup>†</sup>	‡
28	0.26 <sup>†</sup>	0.25 <sup>†</sup>
<b>D8S1179</b>		
8	‡	1.75
9	0.51 <sup>†</sup>	1.00 <sup>†</sup>
10	3.08	8.00
11	4.36	6.25
12	10.51	14.25
13	19.74	34.75
14	33.33	18.75
15	21.03	13.00
16	6.41	2.00
17	1.03 <sup>†</sup>	0.25 <sup>†</sup>
18	‡	‡
19	‡	‡

Allele	African-American (n = 195)	U.S. Caucasian (n = 200)
<b>D21S11</b>		
24	‡	‡
24.2	‡	0.50 <sup>†</sup>
25	‡	‡
26	0.26 <sup>†</sup>	‡
27	5.90	3.75
28	21.80	16.25
28.2	‡	‡
29	20.00	20.75
29.2	‡	0.50 <sup>†</sup>
29.3	0.26 <sup>†</sup>	‡
30	16.15	26.25
30.2	2.56	2.50
31	8.97	5.50
31.2	5.90	10.50
32	0.77 <sup>†</sup>	1.25 <sup>†</sup>
32.2	6.92	7.25
33	0.51 <sup>†</sup>	0.25 <sup>†</sup>
33.1	0.26 <sup>†</sup>	‡
33.2	4.10	4.00
34	0.26 <sup>†</sup>	‡
34.2	‡	0.75 <sup>†</sup>
35	3.59	‡
35.2	‡	‡
36	1.54	‡
37	‡	‡
38	0.26 <sup>†</sup>	‡
<b>D18S51</b>		
7	‡	‡
9	‡	‡
10	0.51 <sup>†</sup>	0.50 <sup>†</sup>
10.2	0.51 <sup>†</sup>	‡
11	1.54	2.00
12	5.64	14.25
13	4.10	15.00
13.2	0.51 <sup>†</sup>	‡
14	6.67	16.75
14.2	0.51 <sup>†</sup>	‡

Allele	African-American (n = 195)	U.S. Caucasian (n = 200)
15	17.69	14.25
16	17.44	14.00
17	16.41	10.50
18	11.03	6.00
19	10.26	3.75
20	4.87	1.50
21	1.03 <sup>†</sup>	1.00 <sup>†</sup>
22	1.03 <sup>†</sup>	0.25 <sup>†</sup>
23	0.26 <sup>†</sup>	0.25 <sup>†</sup>
24	‡	‡
25	‡	‡
26	‡	‡
27	‡	‡
<b>D19S433</b>		
9	‡	‡
10	0.77 <sup>†</sup>	‡
10.2	0.26 <sup>†</sup>	‡
11	7.70	0.25 <sup>†</sup>
11.2	0.26 <sup>†</sup>	‡
12	10.51	6.50
12.2	4.10	‡
13	26.92	26.00
13.2	4.62	2.25
14	21.80	32.75
14.2	5.90	3.25
15	7.44	18.25
15.2	4.87	2.50
16	0.77 <sup>†</sup>	5.25
16.2	3.33	2.00
17	‡	0.50 <sup>†</sup>
17.2	0.51 <sup>†</sup>	0.25 <sup>†</sup>
18.2	0.26 <sup>†</sup>	0.25 <sup>†</sup>
<b>TH01</b>		
4	‡	‡
5	0.26 <sup>†</sup>	‡
6	13.08	25.25
7	40.00	16.00
8	22.82	9.25

Allele	African-American (n = 195)	U.S. Caucasian (n = 200)
9	13.08	13.75
9.3	10.00	35.00
10	0.77 <sup>†</sup>	0.75 <sup>†</sup>
11	‡	‡
13.3	‡	‡
<b>FGA</b>		
16.2	0.26 <sup>†</sup>	‡
17	0.26 <sup>†</sup>	‡
18	0.51 <sup>†</sup>	1.50
19	4.62	6.25
19.2	0.26 <sup>†</sup>	‡
20	4.36	16.25
20.2	‡	0.75 <sup>†</sup>
21	13.33	17.75
21.2	0.26 <sup>†</sup>	‡
22	18.97	16.50
22.2	0.26 <sup>†</sup>	0.50 <sup>†</sup>
23	18.97	14.00
24	16.41	13.25
25	12.31	11.25
26	4.10	1.50
26.2	‡	‡
27	4.10	0.50 <sup>†</sup>
28	0.77 <sup>†</sup>	‡
29	0.26 <sup>†</sup>	‡
30	‡	‡
30.2	‡	‡
31.2	‡	‡
32.2	‡	‡
33.2	‡	‡
34.2	‡	‡
42.2	‡	‡
43.2	‡	‡
44.2	‡	‡
45.2	‡	‡
46.2	‡	‡
47.2	‡	‡
48.2	‡	‡

Allele	African-American (n = 195)	U.S. Caucasian (n = 200)
50.2	‡	‡
51.2	‡	‡

† A minimum allele frequency of 1.3% is suggested by the National Research Council in forensic calculations using either of these African American or U.S. Caucasian databases analyzed using the SGM Plus™ Kit.

‡ Present in the AmpFSTR™ SGM Plus™ Allelic Ladder but not observed in the respective population database samples.

### Analyzing the two databases

Analysis across both databases of 790 total chromosomes (per locus) revealed a total of 11 different D3S1358 alleles, 11 different vWA alleles, 9 different D16S539, 14 different D2S1338 alleles, 10 different D8S1179 alleles, 21 different D21S11 alleles, 17 different D18S51 alleles, 17 D19S433 alleles, 7 different TH01 alleles, and 18 different FGA alleles.

In addition to the alleles that were observed and recorded in the Applied Biosystems databases, other known alleles (listed in [Table 1, “SGM Plus™ Kit loci and alleles” on page 12](#)) have either been published or reported to us by other laboratories.

### Independent allele frequencies

Independence of allelic frequencies within a locus can be expressed by the Hardy-Weinberg (HW) relationship. Approximation of HW expectations in a sample population allows estimation of genotypic frequencies (HW proportions) from observed allelic frequencies using the HW equation (expanded binomial square law) (Hartl and Clark, 1989; Weir, 1996).

Several biostatistical tests were used to survey HW relationships at the SGM Plus™ Kit STR loci in each sample population. Independence was found between alleles within each locus, as  $p$  values  $>0.05$  were obtained from the homozygosity test (Chakraborty *et al.*, 1988; Nei and Roychoudhury, 1974; Nei, 1978), likelihood-ratio test (Edwards *et al.*, 1992; Weir, 1992), and Guo-Thompson exact test (Guo and Thompson, 1992).

Additionally, allele frequency data were analyzed for independence based on the total number of observed distinct homozygous and heterozygous genotype classes (Nei, 1978; Chakraborty *et al.*, 199). Observed values were within two standard errors of expected values for each locus. These sets of data demonstrate that appropriate estimations of SGM Plus™ Kit genotype frequencies are generated from allele frequencies observed in the Life Technologies African-American and U.S. Caucasian databases.

### Random association

Existence of random association (linkage equilibrium) between all 10 STR loci was established through two separate statistical tests. Results of the first test, which considers the observed variance of the number of heterozygous loci (Brown *et al.*, 1980; Budowle *et al.*, 1995) indicate that in both population samples, all SGM Plus™ Kit loci are inherited independently.

Pairwise interclass correlation tests were performed between every possible two-locus combination across the African-American and U.S. Caucasian databases (Karlin *et al.*, 1981). Mendelian behavior between the ten STR loci was observed. SGM Plus™ Kit multilocus genotype frequency estimates may be derived through direct multiplication of each single-locus genotype frequency (the “product rule”) estimated from the Applied Biosystems African-American and U.S. Caucasian databases.

### Low frequency alleles

Some alleles of the SGM Plus™ Kit loci occur at a low frequency (less than five times in either database). For these alleles, a minimum frequency of 0.013 (five divided by 2n, where n equals the number of individuals in the database) was assigned for the SGM Plus™ Kit African-American and U.S. Caucasian databases, as suggested in the 1996 report of the Committee on DNA Forensic Science (National Research Council, 1996). These databases are summarized in [Table 4, “SGM Plus™ Kit allele frequencies” on page 93](#). The minimum reportable genotype frequency at each locus is then  $1.69 \times 10^{-4}$ , giving a minimum combined multilocus genotype frequency of  $1.90 \times 10^{-38}$  for both the African-American and U.S. Caucasian databases.

## Mutation rate

Estimation of spontaneous or induced germline mutation at genetic loci can be achieved by comparing the genotypes of offspring to those of their parents. From such comparisons the number of observed mutations are counted directly.

Genotypes of the ten STR loci amplified by the SGM Plus™ Kit were determined for a total of 146 parent-offspring allelic transfers (meioses) at the Forensic Science Service, Birmingham, England. One length-based STR mutation was observed at the D18S11 locus; mutation was not detected at any of the other nine STR loci. The D18S11 mutation was represented by an increase of one 4-bp repeat unit, a 17 allele was inherited as an 18 (single-step mutation). The maternal/paternal source of this mutation could not be distinguished.

Additional studies (Edwards *et al.*, 1991; Edwards *et al.*, 1992; Weber and Wong, 1993; Hammond *et al.*, 1994; Brinkmann *et al.*, 1995; Chakraborty *et al.*, 1996; Chakraborty *et al.*, 1997; Brinkmann *et al.*, 1998; Momhinweg *et al.*, 1998; Szibor *et al.*, 1998) of direct mutation rate counts produced:

- Larger sample sizes for some of the SGM Plus™ Kit loci.
- Methods for modifications of these mutation rates (to infer mutation rates indirectly for those loci where the rates are not large enough to be measured directly and/or to account for those events undetectable as Mendelian errors).

## Probability of identity

Table 5 shows the Probability of Identity ( $P_I$ ) values of the SGM Plus™ Kit loci individually and combined.

Table 5 Probability of Identity values for the SGM Plus™ Kit STR loci

Locus	African-American	U.S. Caucasian
D3S1358	0.102	0.078
vWA	0.058	0.065
D16S539	0.066	0.103
D2S1338	0.021	0.024
D8S1179	0.075	0.067
D21S11	0.033	0.045
D18S51	0.028	0.030
D19S433	0.039	0.078
TH01	0.102	0.094
FGA	0.035	0.036
Combined	$7.91 \times 10^{-14}$	$2.99 \times 10^{-13}$

The  $P_I$  value is the probability that two individuals selected at random will have an identical SGM Plus™ Kit genotype (Sensabaugh, 1982). The  $P_I$  values for the populations described in this section are then approximately  $1/1.3 \times 10^{13}$  (African-American) and  $1/3.3 \times 10^{12}$  (U.S. Caucasian).

Of 18,915 and 19,900 pairs of SGM Plus™ Kit profiles represented by the African-American and U.S. Caucasian databases, respectively, no 10-locus matches were observed.

Linkage disequilibrium between the SGM Plus™ Kit loci and unique loci included in the AmpFSTR™ Profiler™ PCR Amplification Kit (TPOX, CSF1PO, D5S818, D13S317, and D7S820) was not detected. The combination of these 15 AmpFSTR™ kit loci offers an average probability of identity of approximately  $1/2.2 \times 10^{18}$  (African-American) and  $1/2.1 \times 10^{17}$  (U.S. Caucasian).

## Probability of paternity exclusion

Table 6 shows the Probability of Paternity Exclusion ( $P_E$ ) values of the SGM Plus™ Kit STR loci individually and combined.

Table 6 Probability of paternity exclusion values for the SGM Plus™ Kit loci

Locus	African-American	U.S. Caucasian
D3S1358	0.5260	0.5797
vWA	0.6394	0.6170
D16S539	0.6146	0.5252
D2S1338	0.7818	0.7678

Locus	African-American	U.S. Caucasian
D8S1179	0.5930	0.6128
D21S11	0.7281	0.6835
D18S51	0.7518	0.7414
D19S433	0.7031	0.5865
TH01	0.5250	0.5418
FGA	0.7202	0.7173
Combined	0.99998	0.99997

The  $P_E$  value is the probability, averaged over all possible mother-child pairs, that a random alleged father will be excluded from paternity after DNA typing of the SGM Plus™ Kit STR loci (Chakraborty and Stivers, 1996).

The  $P_E$  value offered by the combination of the 10 AmpF $\phi$ STR™ kit loci (D3S1358, vWA, D16S539, D2S1338, D8S1179, D21S11, D18S51, D19S433, TH01, FGA) amplified by the AmpF $\phi$ STR SGM Plus and Profiler kits is approximately 0.9999997 for the African American database and approximately 0.9999992 for the U.S. Caucasian database.

## Section 5.2 Performance Validation After Buffer and Enzyme Component Replacement

### Overview

As part of an ongoing program to exercise greater control over raw materials used in the AmpF $\ell$ STR<sup>™</sup> PCR Amplification Kits, manufacturing of the AmpliTaq Gold<sup>™</sup> enzyme and 10X PCR Buffer II (Tris-KCl buffer) components is transitioning from Roche Molecular Systems to Life Technologies. Manufacturing of both components by Life Technologies will be conducted according to the same specifications used previously by Roche. The in-house components are established raw materials in our next generation kits (for example, the NGM<sup>™</sup>, NGM SElect<sup>™</sup> and Identifiler<sup>™</sup> Plus Kits).

### Experiments

We performed studies to compare the performance of the SGM Plus<sup>™</sup> Kit containing the in-house components (updated kit) with the performance of the original kit, focussing on studies most relevant to forensic DNA testing (see SWGDAM Guidelines effective January 1, 2011). These studies, while not exhaustive, are in our opinion appropriate for a manufacturer.

Our studies compared the performance of two Roche-manufactured enzyme and buffer lots (Control mixes) with three new lots of buffer and two new lots of enzyme manufactured by Life Technologies (Test mixes). Studies were performed using Test mixes containing both the enzyme and buffer manufactured by Life Technologies.

Test Material	Control A mix	Control B mix	Test A mix	Test B mix	Test C mix
Buffer	Control Buffer Lot 1	Control Buffer Lot 2	Test Buffer Lot 1	Test Buffer Lot 2	Test Buffer Lot 3
Enzyme	Control Enzyme Lot 1	Control Enzyme Lot 2	Test Enzyme Lot 1	Control Enzyme Lot 2	Test Enzyme Lot 2

Each of the five mixes listed above were used to conduct reproducibility, sensitivity, and inhibition studies. All amplifications were performed using a GeneAmp<sup>™</sup> PCR System 9700 with either silver or gold-plated silver block using the recommended amplification conditions and cycle number for the SGM Plus<sup>™</sup> Kit. All data was run on an Applied Biosystems 3130xl Genetic Analyzer running Data Collection Software v3.0 and analyzed using GeneMapper<sup>™</sup> ID-X Software. Subsequent data analysis was performed using Minitab<sup>™</sup> Statistical Software.

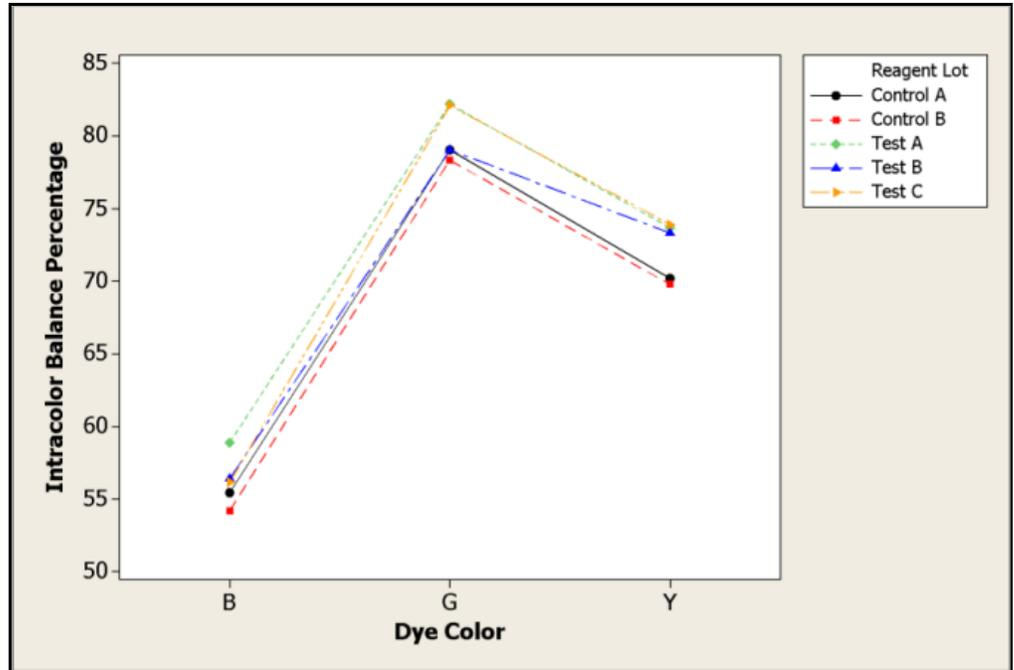
## Reproducibility study

For the reproducibility study, 12 replicates of control DNA 007 at 2 ng input and three negative control replicates were amplified. The results were evaluated for intracolor balance, stutter percentage, and the presence, signal intensity, and location of artifacts.

### Intracolor balance

No significant difference (<10% increase or decrease) in the level of intracolor balance was observed between the Test and Control mixes (Figure 17).

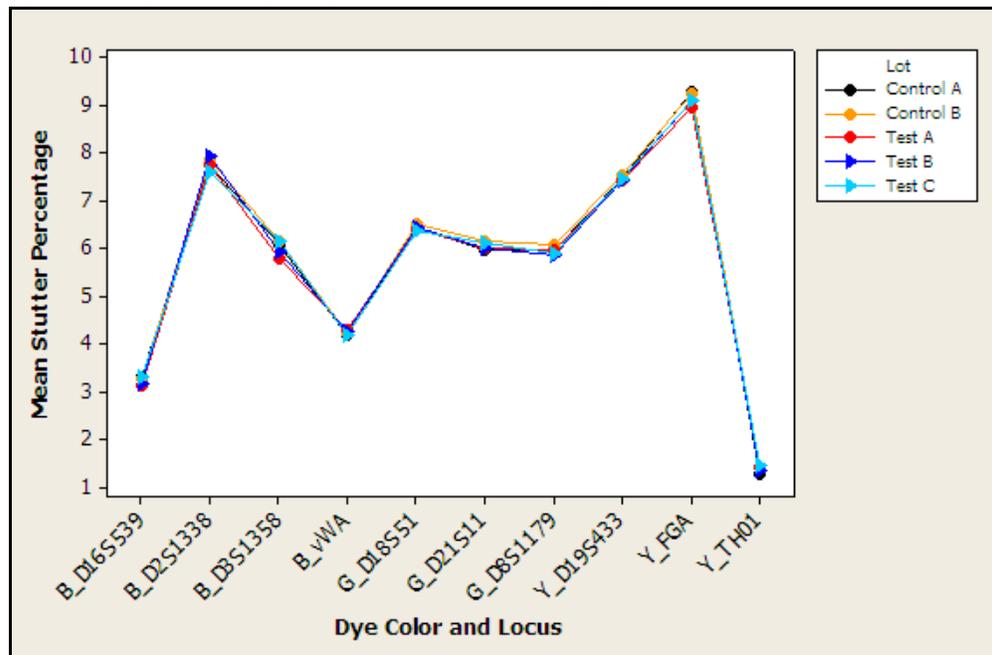
Figure 17 Reproducibility study: intracolor balance



### Stutter percentages

Stutter percentage results for each marker were comparable across all Test and Control mixes (Figure 18).

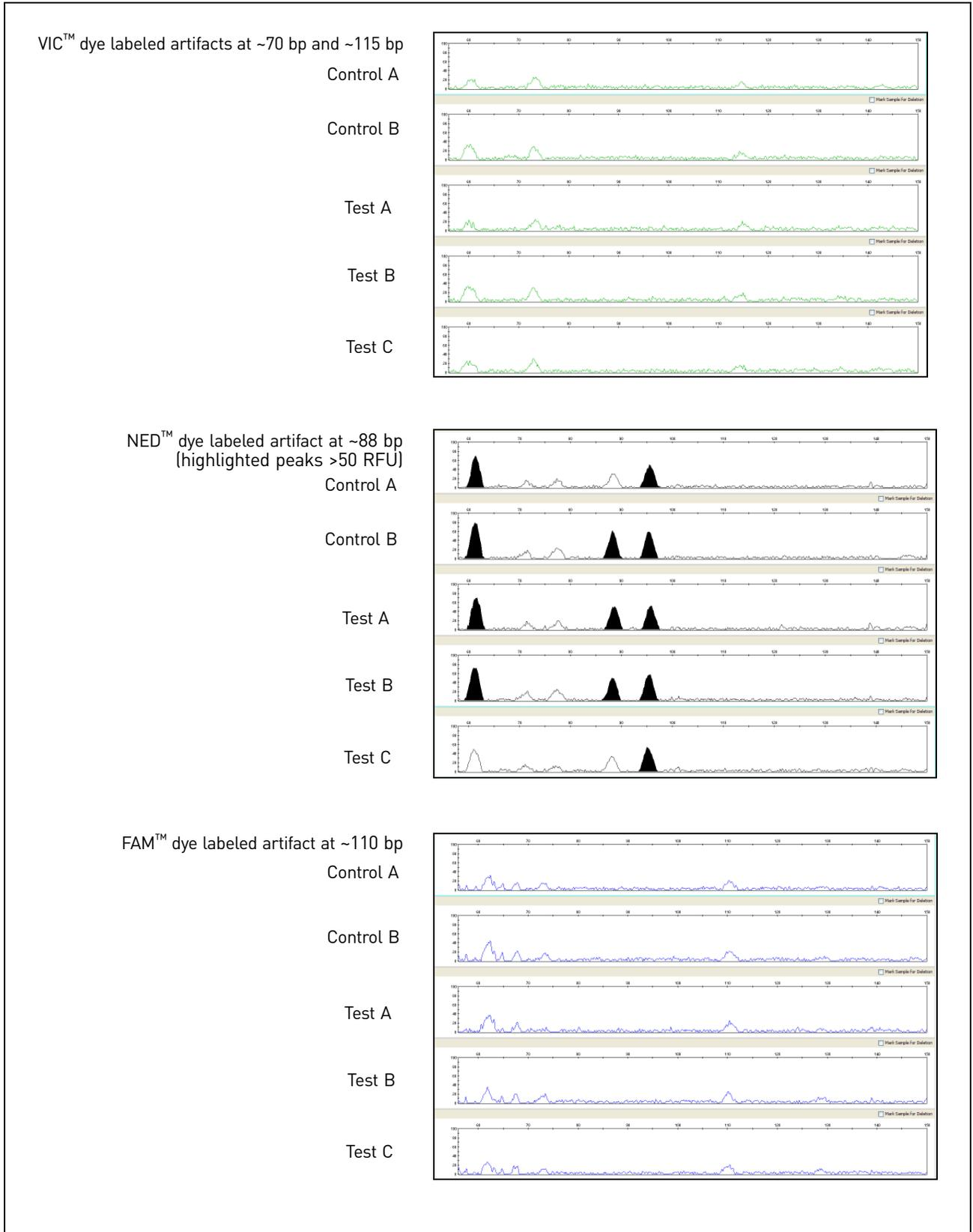
Figure 18 Reproducibility study: mean stutter percentage



## Artifacts

Known artifacts observed showed the same morphology, signal intensity, and location in all Test and Control mixes and do not exceed 50 RFU, except for the NED™ dye artifacts highlighted in Figure 19. No new artifacts were observed in the Test mixes.

Figure 19 Reproducibility study: known artifacts (Y-scale 100 RFU)



Performance Validation: Buffer and Enzyme Replacement

## Sensitivity study

For the sensitivity study, dilution series of three genomic DNA samples were amplified: 2 ng (three replicates), 1 ng, 0.5 ng, and 0.25 ng (four replicates each). The results were evaluated for mean peak height, degree of linearity between input DNA concentration and peak height, level of allelic dropout at 250 pg, and genotype concordance.

### Mean peak height

Overall mean peak height observations were consistent between all Test and Control mixes (Figure 20) demonstrating equivalent performance (Figure 21). Sample IBB286 produced a wider spread between data points. However, this spread did not correlate to a difference in Test versus Control mixes and is likely due to variability in sample set up.

Figure 20 Sensitivity study: mean peak heights for three genomic DNA samples

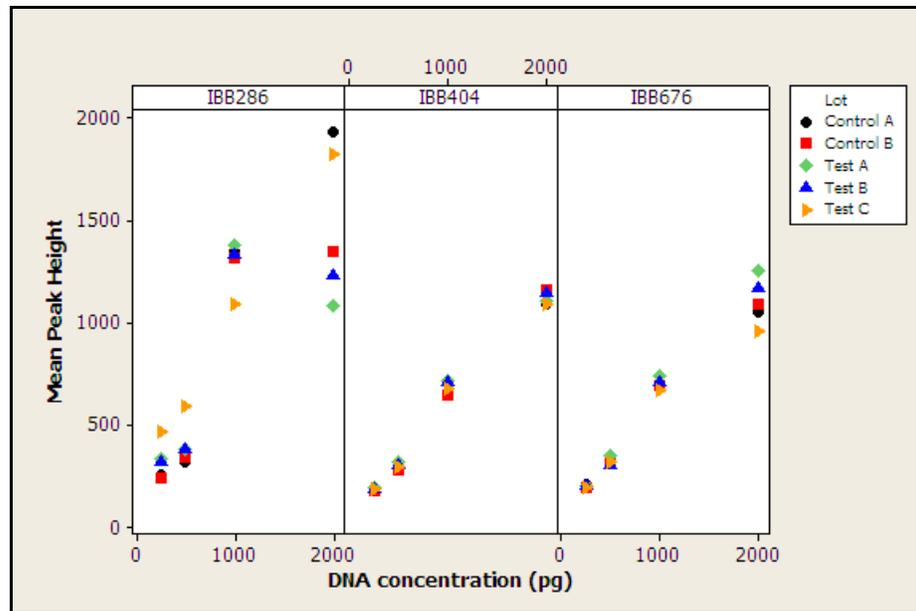
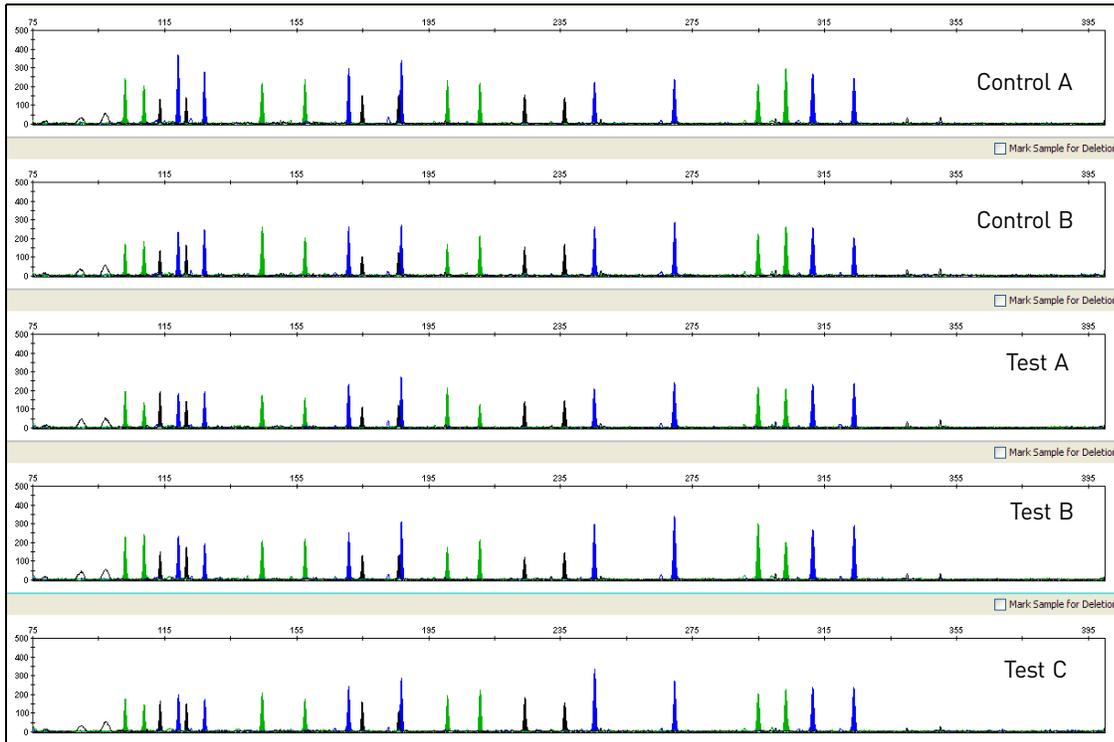


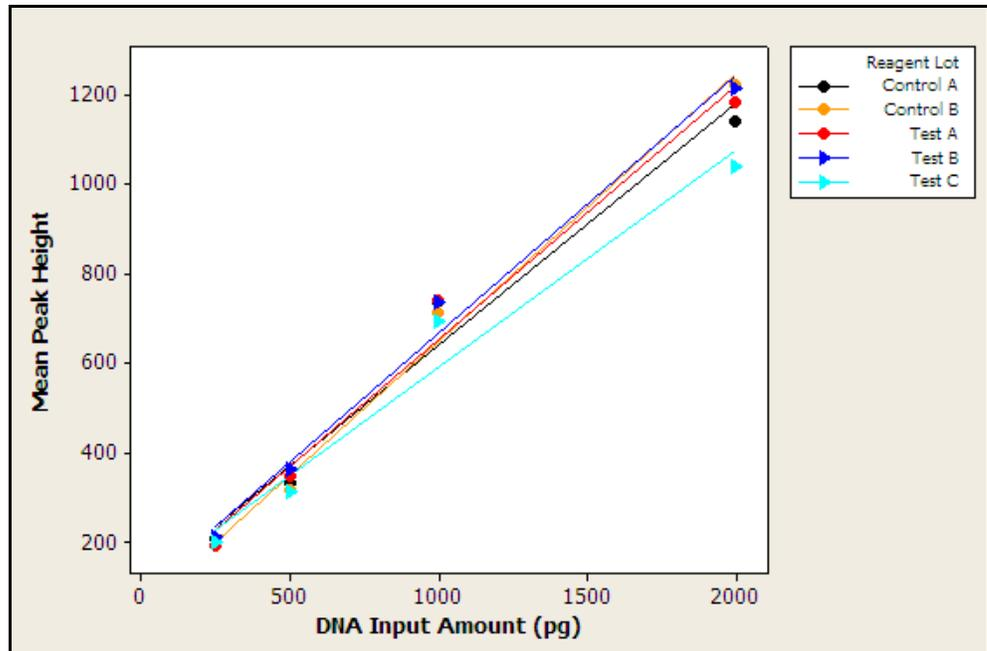
Figure 21 Sensitivity study: representative electropherograms for Sample 2 amplified using 250 pg input DNA (Y-scale 500 RFU)



**DNA concentration and peak height**

The calculated slope and  $R^2$  values for each of the plotted curves are equivalent, showing comparable relationships between peak height and DNA input amount for the Test and Control mixes (Figure 22).

Figure 22 Sensitivity study: linear regression plot of combined mean peak height for three genomic DNA samples



**Allelic dropout**

No allelic dropout was observed for any of the Test or Control mixes.

**Genotype concordance**

Genotypes for Test and Control mixes were 100% concordant (Table 7).

**Table 7** Sensitivity study: genotype concordance

DNA Input Amount	Reagent Mix	Genotype Concordance
250 pg	Test A	100%
	Test B	100%
	Test C	100%
	Control A	100%
	Control B	100%
500 pg	Test A	100%
	Test B	100%
	Test C	100%
	Control A	100%
	Control B	100%
1 ng	Test A	100%
	Test B	100%
	Test C	100%
	Control A	100%
	Control B	100%
2 ng	Test A	100%
	Test B	100%
	Test C	100%
	Control A	100%
	Control B	100%

**Inhibition study**

An inhibition series of 2 ng control DNA 007 consisting of uninhibited control, humic acid at a final concentration of 15.2 ng/ $\mu$ L, and hematin at a final concentration of 40  $\mu$ M in replicates of five were amplified using each of Test and Control mixes. The amount of each inhibitor tested was titrated to cause an approximate 50% reduction in overall peak height of the samples. Results were evaluated for mean peak height, minimum peak height, intracolor balance, and levels of allelic dropout.

**Mean peak height, minimum mean peak height, and intracolor balance**

No significant difference in mean peak height, mean minimum peak height and intracolor balance was observed for any Test or Control mixes tested on uninhibited Control DNA 007 and Control DNA 007 inhibited with hematin or humic acid. (Figure 23, 24, and 25).

Figure 23 Inhibition study: mean peak height

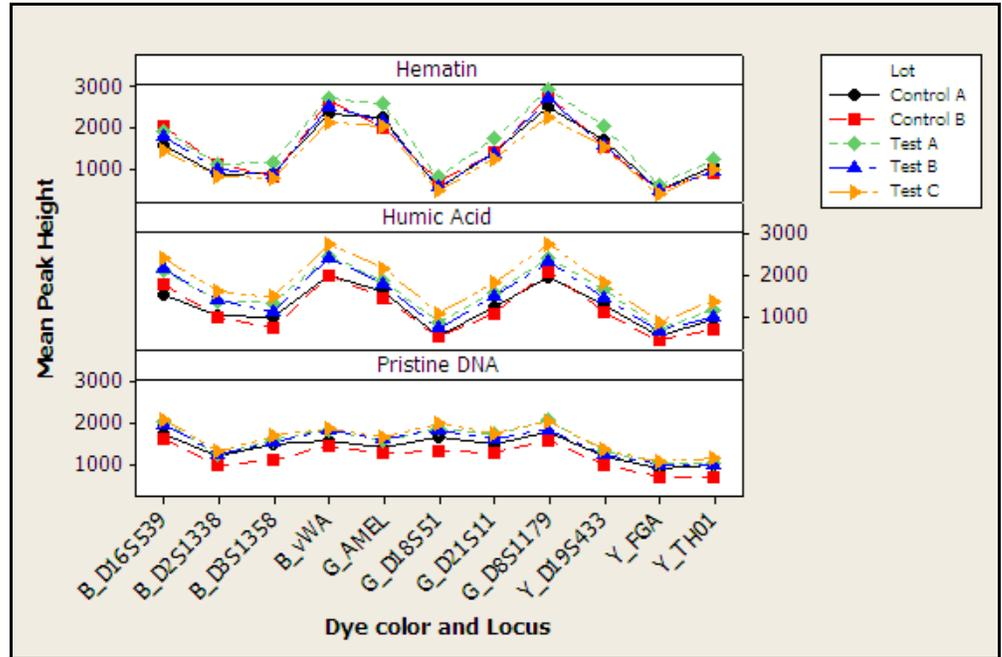


Figure 24 Inhibition study: minimum peak height

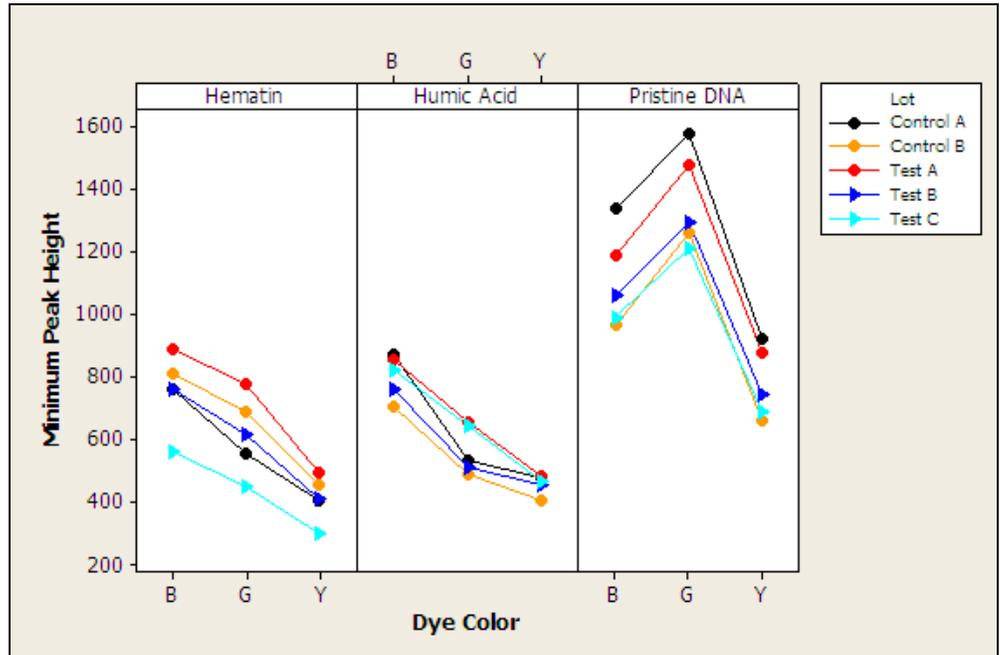
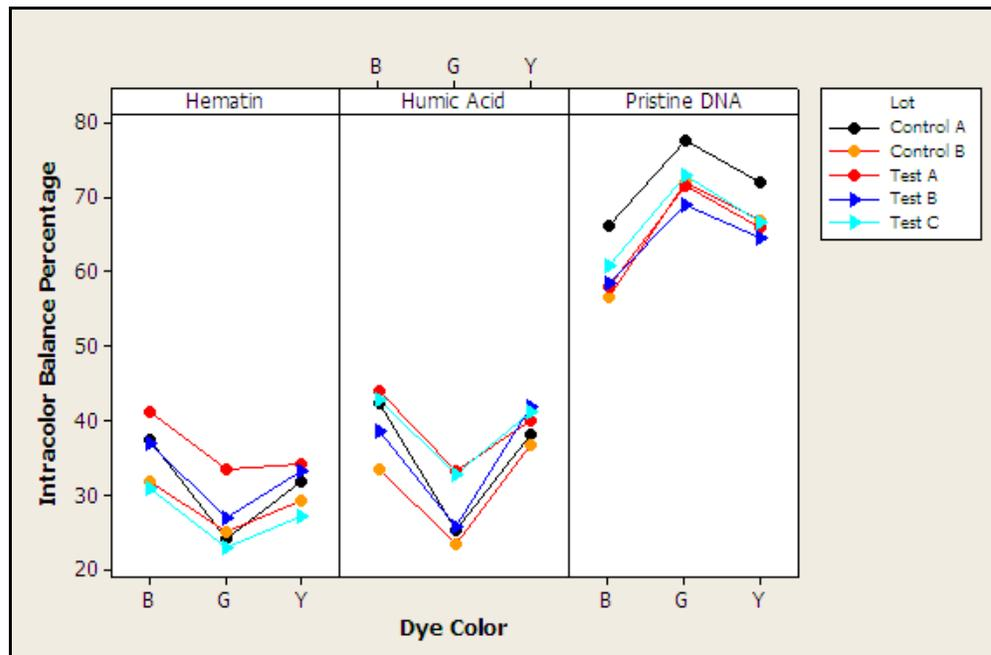
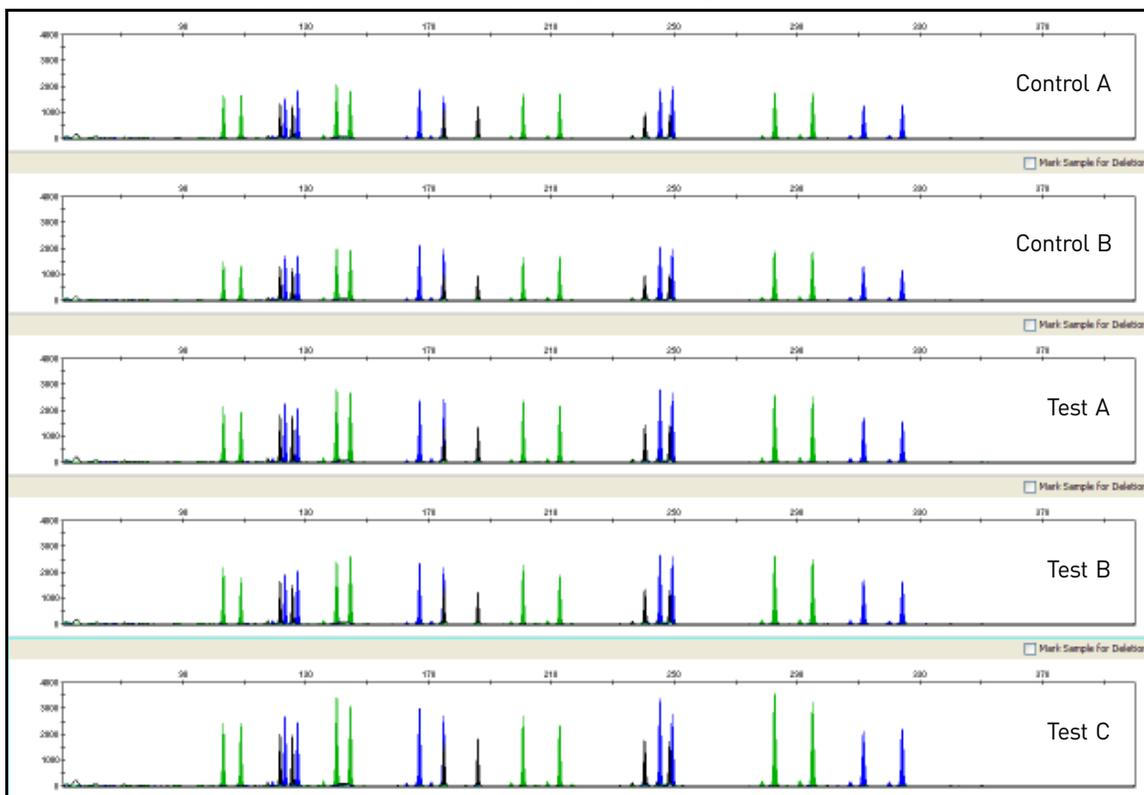


Figure 25 Inhibition study: intracolor balance

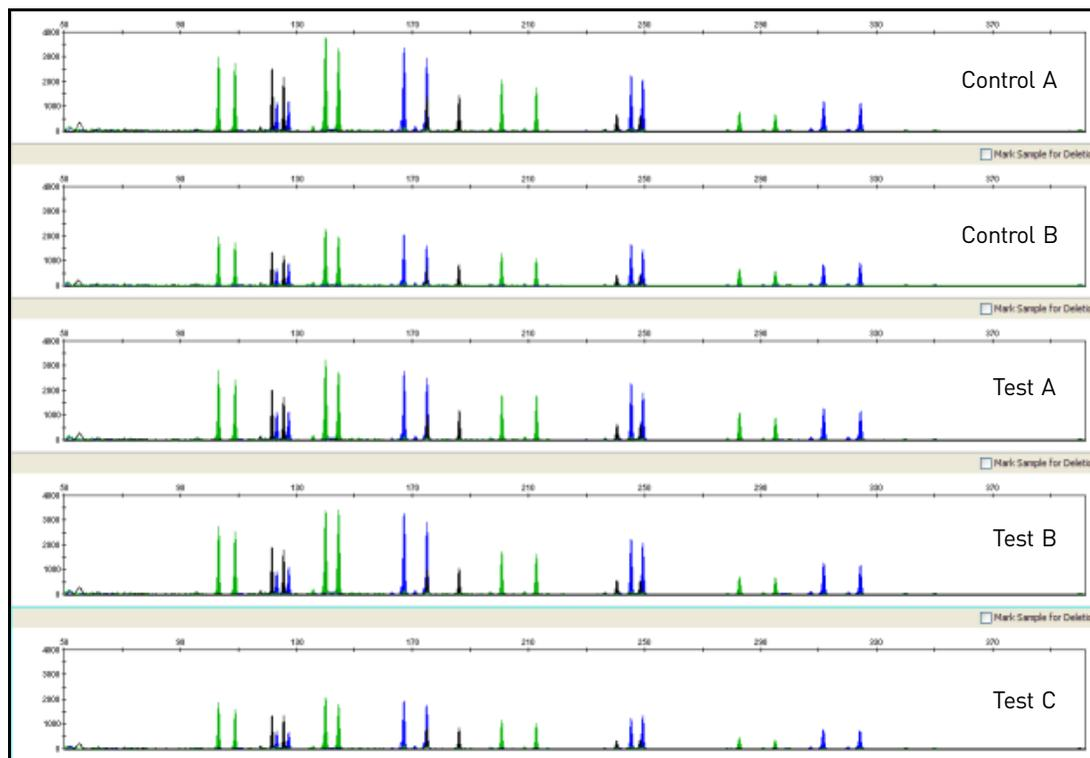


Representative electropherograms from the inhibition study are shown in [Figure 26](#), [Figure 27](#), and [Figure 28](#).

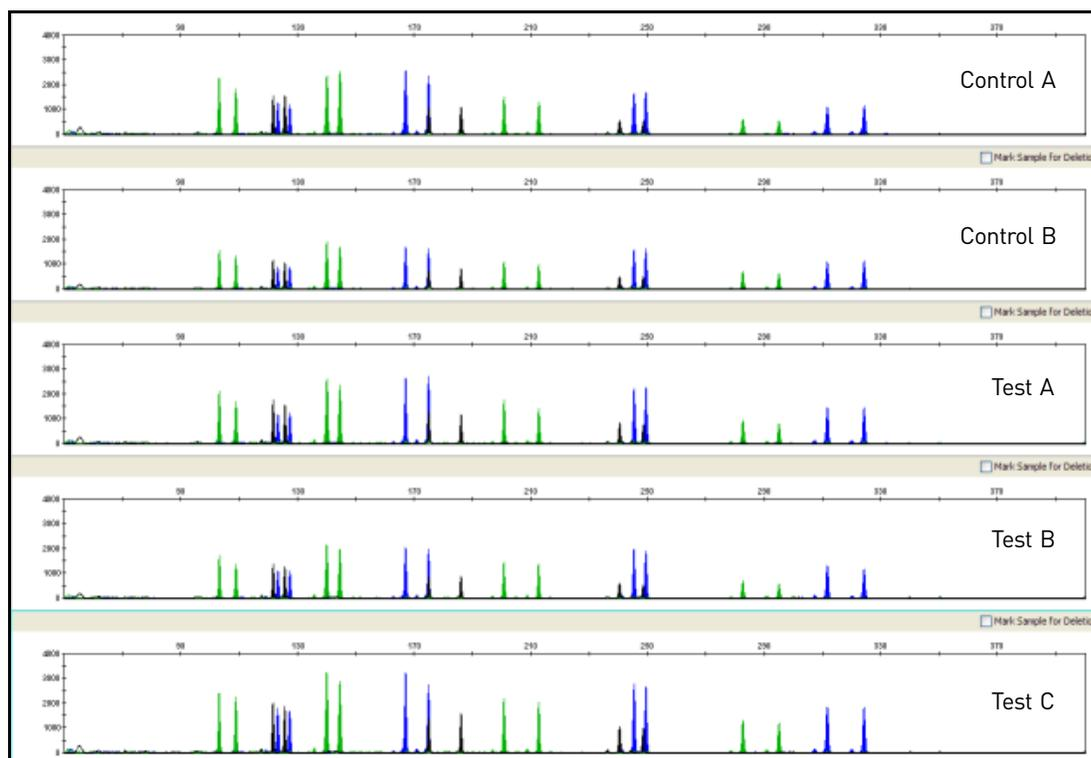
Figure 26 Inhibition study: representative electropherograms using uninhibited Control DNA 007 (Y-scale 4000 RFU)



**Figure 27** Inhibition study: representative electropherograms using Control DNA 007 inhibited with 40  $\mu$ M Hematin (Y-scale 4000 RFU)



**Figure 28** Inhibition study: representative electropherograms using Control DNA 007 inhibited with 15.25 ng/ $\mu$ L Humic Acid (Y-scale 4000 RFU)



**Allelic dropout**

No allelic dropout events were seen for any Test or Control mixes tested on uninhibited Control DNA 007 and Control DNA 007 inhibited with hematin or humic acid.

**Conclusions**

Laboratories can expect to obtain equivalent quality profiles across a wide range of forensic samples when using the SGM Plus™ Kit containing the AmpliTaq Gold™ enzyme and 10X PCR Buffer II manufactured by Life Technologies as compared to the original SGM Plus™ Kit containing AmpliTaq Gold™ enzyme and 10X PCR Buffer II manufactured by Roche Molecular Systems.



# Troubleshooting

Follow the actions recommended in this appendix to troubleshoot problems that occur during analysis.

**Table 8** Troubleshooting

Observation	Possible causes	Recommended actions
Faint or no signal from both the AmpF $\mathcal{L}$ STR™ Control DNA 007 and the DNA test samples at all loci	Incorrect volume or absence of AmpF $\mathcal{L}$ STR™ PCR Reaction Mix, SGM Plus™ Primer Set, or AmpliTaq Gold™ DNA Polymerase	Repeat amplification.
	No activation of AmpliTaq Gold™ DNA Polymerase	Repeat amplification, making sure to hold reactions initially at 95°C for 11 minutes.
	Master Mix not vortexed thoroughly before aliquoting	Vortex the Master Mix thoroughly.
	SGM Plus™ Primer Set exposed to too much light	Store the Primer Set protected from light.
	GeneAmp™ PCR System malfunction	Refer to the thermal cycler user's manual and check instrument calibration.
	Use of incorrect thermal cycling parameters	Check the protocol for correct thermal cycling parameters.
	Tubes not seated tightly in the thermal cycler during amplification	Push reaction tubes firmly into contact with block after first cycle. Repeat test.
	Wrong PCR reaction tube	Use Applied Biosystems MicroAmp™ Reaction Tubes with Caps or a MicroAmp™ optical 96-well plate.
	MicroAmp™ Base used with tray/retainer set and tubes in GeneAmp™ 9700	Remove MicroAmp™ Base from tray/retainer set and repeat test.
	Insufficient PCR product electrokinetically injected	Prepare PCR product as described in <a href="#">Chapter 3, "Electrophoresis"</a> on page 25.
Degraded formamide	Check the storage of formamide; do not thaw and refreeze multiple times. Try Hi-Di™ Formamide.	

Observation	Possible causes	Recommended actions
Positive signal from AmpF $\lambda$ STR™ Control DNA 007 but partial or no signal from DNA test samples	Quantity of test DNA sample is below assay sensitivity	Quantitate DNA and add 1.0–2.5 ng of DNA. Repeat test.
	Test sample contains high concentration of PCR inhibitor (for example, heme compounds, certain dyes)	Quantitate DNA and add minimum necessary volume. Repeat test.
		Wash the sample in a Centricon™-100 centrifugal filter unit. Repeat test.
	Test sample DNA is severely degraded	If possible, evaluate the quality of DNA sample by running an agarose gel. If DNA is degraded, reamplify with an increased amount of DNA or use the AmpF $\lambda$ STR™ MiniFiler™ Kit.
	Dilution of test sample DNA in water or wrong buffer (for example, TE formula with incorrect EDTA concentration)	Redilute DNA using low-TE Buffer (with 0.1 mM EDTA).
More than one allele present at a locus	Presence of exogenous DNA	Use appropriate techniques to avoid introducing foreign DNA during laboratory handling.
	Amplification of stutter product	See <a href="#">"Stutter products"</a> on page 74.
	Mixed sample	
Some but not all loci visible on electropherogram of DNA test samples	Test-sample DNA is severely degraded	If possible, evaluate the quality of DNA sample by running an agarose gel. If DNA is degraded, reamplify with an increased amount of DNA or use the AmpF $\lambda$ STR™ MiniFiler™ Kit.
	Test sample contains high concentrations of a PCR inhibitor (for example, heme compounds, certain dyes)	Quantitate DNA and add minimum necessary volume. Repeat test.
		Wash the sample in a Centricon™-100 centrifugal filter unit. Repeat test.
Poor peak height balance	Incorrect thermal cycler parameters	Check the protocol for correct thermal cycler parameters.
	GeneAmp™ PCR System 9700 with Aluminum 96-Well block or third-party thermal cyclers	Use Applied Biosystems GeneAmp™ PCR System 9700 with silver or gold-plated silver blocks only, or the Veriti™ 96-Well Thermal Cycler.



# PCR Work Areas

- Work area setup and lab design . . . . . 115
- PCR setup work area . . . . . 115
- Amplified DNA work area . . . . . 116

## Work area setup and lab design

Many resources are available for the appropriate design of a PCR laboratory. If you are using the AmpF $\Phi$ STR™ SGM Plus™ PCR Amplification Kit for:

- Forensic DNA testing, refer to “Forensic Laboratories: Handbook for Facility Planning, Design, Construction and Moving,” National Institute of Justice, 1998
- Parentage DNA testing, refer to the “Guidance for Standards for Parentage Relationship Testing Laboratories,” American Association of Blood Banks, 7th edition, 2004

The sensitivity of the AmpF $\Phi$ STR™ Kits (and other PCR-based tests) enables amplification of minute quantities of DNA, necessitating precautions to avoid contamination of samples yet to be amplified (Kwok and Higuchi, 1989).

Also take care while handling and processing samples to prevent contamination by human DNA. Wear gloves at all times and change them frequently. Close sample tubes when not in use. Limit aerosol dispersal by handling sample tubes and reagents carefully.

**Note:** We do not intend these references for laboratory design to constitute all precautions and care necessary for using PCR technology.

## PCR setup work area

---

**IMPORTANT!** These items should never leave the PCR Setup Work Area.

---

- Calculator
- Gloves, disposable
- Marker pen, permanent
- Microcentrifuge
- Microcentrifuge tubes, 1.5-mL, or 2.0-mL, or other appropriate clean tube (for Master Mix preparation)
- Microcentrifuge tube rack
- Pipette tips, sterile, disposable hydrophobic filter-plugged
- Pipettors

- Tube decapper, autoclavable
- Vortex

## Amplified DNA work area

---

**IMPORTANT!** Place the thermal cyclers in the Amplified DNA Work Area.

---

You can use the following systems:

- GeneAmp™ PCR System 9700 with the Silver 96-Well Block
- GeneAmp™ PCR System 9700 with the Gold-plated Silver 96-Well Block

---

**IMPORTANT!** The SGM Plus™ Kit is not validated for use with the GeneAmp™ PCR System 9700 with the Aluminium 96-Well Block. Use of this thermal cycling platform may adversely affect performance of the kit.

---

- Veriti™ 96-Well Thermal Cycler



# Ordering Information

## Equipment and materials not included

Table 9 and Table 10 list required and optional equipment and materials not supplied with the SGM Plus™ Kit. Unless otherwise noted, many of the items are available from major laboratory suppliers (MLS).

Table 9 Equipment

Equipment	Source
Applied Biosystems 3100/3100- <i>Avant</i> Genetic Analyzer	Contact your local Life Technologies sales representative
Applied Biosystems 3130/3130 <i>xl</i> Genetic Analyzer	
Applied Biosystems 3500/3500 <i>xL</i> Genetic Analyzer for Human Identification	
Applied Biosystems 310 Genetic Analyzer	
GeneAmp™ PCR System 9700 with the Silver 96-Well Block	N8050001
GeneAmp™ PCR System 9700 with the gold-plated silver 96-well block	4314878
Veriti™ 96-Well Thermal Cycler	4375786
Silver 96-well sample block	N8050251
Gold-plated silver 96-well sample block	4314443
Tabletop centrifuge with 96-well plate adapters (optional)	MLS

Table 10 User-supplied materials

Item†	Source
AmpF <del>L</del> STR™ SGM Plus™ PCR Amplification Kit	4307133
<b>3100 Analyzer materials</b>	
96-well plate septa	4315933
Reservoir septa	4315932
3100/3130 <i>xl</i> Genetic Analyzer capillary array, 36-cm	4315931
POP-4™ polymer for 3100/3100- <i>Avant</i> Genetic Analyzers	4316355
3100/3100- <i>Avant</i> Genetic Analyzer Autosampler Plate Kit, 96-well	4316471
GeneScan™ 500 ROX™ Size Standard	401734
Running Buffer, 10X	402824
Hi-Di™ Formamide	4311320
DS-32 Matrix Standard Kit (Dye Set F)	4345831
MicroAmp™ Optical 96-well reaction plate	N8010560



Item <sup>†</sup>	Source
250-µL glass syringe (array-fill syringe)	4304470
5.0-mL glass syringe (polymer-reserve syringe)	628-3731
For a complete list of parts and accessories for the 3100/3100- <i>Avant</i> instrument, refer to Appendix B of the <i>3100 Genetic Analyzer and 3100-<i>Avant</i> Genetic Analyzer User Reference Guide</i> (Pub. no. 4335393).	
<b>3130xl Analyzer materials</b>	
96-well plate septa	4315933
Reservoir septa	4315932
3100/3130xl Genetic Analyzer capillary array, 36-cm	4315931
POP-4™ polymer for 3130/3130xl Genetic Analyzers	4352755
3100/3100- <i>Avant</i> Genetic Analyzer Autosampler Plate Kit, 96-well	4316471
GeneScan™ 500 ROX™ Size Standard	401734
Running Buffer, 10X	402824
DS-32 Matrix Standard Kit (Dye Set F)	4345831
MicroAmp™ Optical 96-well reaction plate	N8010560
Hi-Di™ Formamide	4311320
For a complete list of parts and accessories for the 3130/3130xl instrument, refer to Appendix A of the <i>Applied Biosystems 3130/3130xl Genetic Analyzers Maintenance, Troubleshooting, and Reference Guide</i> (Pub. no. 4352716).	
<b>3500/3500xL Analyzer materials</b>	
Anode buffer container (ABC)	4393927
Cathode buffer container (CBC)	4408256
POP-4™ polymer (960 samples) for 3500/3500xL Genetic Analyzers	4393710
POP-4™ polymer (384 samples) for 3500/3500xL Genetic Analyzers	4393715
Conditioning reagent	4393718
8-Capillary array, 36 cm for 3500 Genetic Analyzers	4404683
24-Capillary array, 36 cm for 3500xL Genetic Analyzers	4404687
96-well retainer & base set (Standard) 3500/3500xL Genetic Analyzers	4410228
8-Tube retainer & base set (Standard) for 3500/3500xL Genetic Analyzers	4410231
8-Strip Septa for 3500/3500xL Genetic Analyzers	4410701
96-Well Septa for 3500/3500xL Genetic Analyzers	4412614
Septa Cathode Buffer Container, 3500 series	4410715
GeneScan™ 500 ROX™ Size Standard	401734
For a complete list of parts and accessories for the 3500/3500xL instrument, refer to the <i>Applied Biosystems 3500/3500xL Genetic Analyzer User Guide</i> (Pub. no. 4401661)	
<b>310 Analyzer materials</b>	
310 DNA Analyzer capillary array, 47-cm	402839
0.5 mL sample tray	5572
96-well tray adaptor (for 9700 thermal cycler trays)	4305051
GeneScan™ 500 ROX™ Size Standard	401734

Item <sup>†</sup>	Source
Running Buffer, 10X	4335643
Genetic analyzer septa retainer clips for 96-tube sample tray	402866
Genetic analysis sample tubes (0.5-mL)	401957
Septa for 0.5-mL sample tubes	401956
DS-32 Matrix Standard Kit (Dye Set F) for the 310 Genetic Analyzer	4312131
MicroAmp™ 8-tube strip, 0.2-mL	N8010580
MicroAmp™ 96-well base (holds 0.2-mL reaction tubes)	N8010531
MicroAmp™ 96-well full plate cover	N8010550
MicroAmp™ 96-well tray/retainer set	403081
POP-4™ polymer for the 310 Genetic Analyzer	402838
For a complete list of parts and accessories for the 310 instrument, refer to Appendix B of the <i>310 Genetic Analyzer User Guide</i> (Pub. no. 4317588).	
PCR Amplification	
MicroAmp™ 96-well tray	N8010541
MicroAmp™ reaction tube with cap, 0.2-mL	N8010540
MicroAmp™ 8-tube strip, 0.2-mL	N8010580
MicroAmp™ 8-cap strip	N8010535
MicroAmp™ 96-well tray/retainer set	403081
MicroAmp™ 96-well base	N8010531
MicroAmp™ clear adhesive film	4306311
MicroAmp™ optical adhesive film	4311971
MicroAmp™ optical 96-well reaction plate	N8010560
Other user-supplied materials	
Hi-Di™ Formamide, 25-mL	4311320
Aerosol resistant pipette tips	MLS
Microcentrifuge tubes	MLS
Pipettors	MLS
Tape, labeling	MLS
Tube, 50-mL Falcon	MLS
Tube decapper, autoclavable	MLS
Deionized water, PCR grade	MLS
Tris-HCL, pH 8.0	MLS
EDTA, 0.5 M	MLS
Vortex	MLS

† For the Safety Data Sheet (SDS) of any chemical not distributed by Life Technologies, contact the chemical manufacturer. Before handling any chemicals, refer to the SDS provided by the manufacturer, and observe all relevant precautions.



**Appendix C** Ordering Information  
*Equipment and materials not included*



**WARNING! GENERAL SAFETY.** Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
  - Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, etc). To obtain SDSs, see the “Documentation and Support” section in this document.
- 



## Chemical safety



**WARNING! GENERAL CHEMICAL HANDLING.** To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below, and consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the “Documentation and Support” section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

### Specific chemical handling

CAS	Chemical	Phrase
26628-22-8	Sodium Azide	Sodium azide may react with lead and copper plumbing to form highly explosive metal azides.



## Biological hazard safety



**WARNING! Potential Biohazard.** Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.

---



**WARNING! BIOHAZARD.** Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

In the U.S.:

- U.S. Department of Health and Human Services guidelines published in Biosafety in Microbiological and Biomedical Laboratories found at: [www.cdc.gov/biosafety](http://www.cdc.gov/biosafety)
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030), found at: [www.access.gpo.gov/nara/cfr/waisidx\\_01/29cfr1910a\\_01.html](http://www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html)
- Your company's/institution's Biosafety Program protocols for working with handling potentially infectious materials.
- Additional information about biohazard guidelines is available at: [www.cdc.gov](http://www.cdc.gov)

In the EU:

Check local guidelines and legislation on biohazard and biosafety precaution and refer to the best practices published in the World Health Organization (WHO) Laboratory Biosafety Manual, third edition, found at: [www.who.int/csr/resources/publications/biosafety/WHO\\_CDS\\_CSR\\_LYO\\_2004\\_11/en/](http://www.who.int/csr/resources/publications/biosafety/WHO_CDS_CSR_LYO_2004_11/en/)

---





**Appendix D Safety**  
*Biological hazard safety*



# Bibliography

- Akane, A., Matsubara, K., Nakamura, H., Takahashi, S., and Kimura, K. 1994. Identification of the heme compound copurified with deoxyribonucleic acid (DNA) from bloodstains, a major inhibitor of polymerase chain reaction (PCR) amplification. *J. Forensic Sci.* 39:362–372.
- Barber, M.D., Piercy, R.C., Andersen, J.F., and Parkin, B.H. 1995. Structural variation of novel alleles at the Hum vWA and Hum FES/FPS short tandem repeat loci. *Intl. J. Legal Med.* 108:31–35.
- Andersen, J.F., Greenhalgh M.J., Butler, H.R., Kilpatrick, S.R., Piercy, R.C., Way, K.A., Myhill, H.S., Wright, J.C., Hallett, R., and Parkin, B.H. 1996. Further validation of a multiplex STR system for use in routine forensic identity testing. *Forensic Sci. Int.* 78:47–64.
- Bär, W., Brinkmann, B., Budowle, B., Carracedo, A., Gill, P., Lincoln, P., Mayr, W., and Olaisen, B. 1997. DNA recommendations. Further report of the DNA Commission of the ISFH regarding the use of short tandem repeat systems. *Intl. J. Legal Med.* 110:175–176.
- Barber, M.D., Piercy, R.C., Andersen, J.F., and Parkin, B.H. 1995. Structural variation of novel alleles at the Hum vWA and Hum FES/FPS short tandem repeat loci. *Intl. J. Legal Med.* 108:31–35.
- Barber, M.D., McKeown, B.J., and Parkin, B.H. 1996. Structural variation in the alleles of a short tandem repeat system at the human alpha fibrinogen locus. *Intl. J. Legal Med.* 108:180–185.
- Barber, M.D. and Parkin, B.H. 1996. Sequence analysis and allelic designation of the two short tandem repeat loci D18S51 and D8S1179. *Intl. J. Legal Med.* 109:62–65.
- Begovich A.B., McClure G.R., Suraj V.C., Helmuth R.C., Fildes N., Bugawan T.L., Erlich H.A., and Klitz W. 1992. Polymorphism, recombination, and linkage disequilibrium within the HLA class II region. *J. Immunol.* 148:249–58.
- Brinkmann, B., Moller, A. and Wiegand, P. 1995. Structure of new mutations in 2 STR systems. *Intl. J. Legal Med.* 107:201–203.
- Brinkmann, B., Junge, A., Meyer, E., and Wiegand, P. 1998. Population genetic diversity in relation to microsatellite heterogeneity. *Hum. Mutat.* 11:135-144.
- Brinkmann, B., Klitschar, M., Neuhuber, F., Huhne, J. and Rolf, B. 1998. Mutation rate in human microsatellites: Influence of the structure and length of the tandem repeat. *Am. J. Hum. Genet.* 62:1408–1415.
- Brown, A.H.D., Feldman, M.W., and Nevo, E. 1980. Multilocus structure of natural populations of *Hordeum spontaneum*. *Genetics* 96:523–536.
- Budowle, B. et al. 1995. D1S80 population data in African-Americans, Caucasians, Southeastern Hispanics, Southwestern Hispanics, and Orientals. *J. Forensic Sci.* 40:38–44.

- Buel, E., Wang, G., and Schwartz, M. 1995. PCR amplification of animal DNA with human X-Y amelogenin primers used in gender determination. *J. Forensic Sci.* 40:641–644.
- Chakraborty, R., Stivers, D., and Zhong, Y. 1996. Estimation of mutation rates from parentage exclusion data: applications to STR and VNTR loci. *Mutat. Res.* 354:41–48.
- Chakraborty, R., Smouse, P.E., and Neel, J.V. 1988. Population amalgamation and genetic variation: observations on artificially agglomerated tribal populations of Central and South America. *Am. J. Hum. Genet.* 43:709–725.
- Chakraborty, R., Fornage, M., Guegue, R., and Boerwinkle, E. 1991. Population genetics of hypervariable loci: analysis of PCR based VNTR polymorphism within a population. In: Burke, T., Doif, G., Jeffreys, A.J., and Wolff, R., eds. *DNA Fingerprinting: Approaches and Applications*. Birkhauser Verlag, Berlin, pp. 127–143.
- Chakraborty, R. and Stivers, D.N. 1996. Paternity exclusion by DNA markers: effects of paternal mutations. *J. Forensic Sci.* 41:671–677.
- Chakraborty, R., Kimmel, M., Stivers, D., Davison, L., and Deka, R. 1997. Relative mutation rates at di-, tri-, and tetranucleotide microsatellite loci. *Proc. Natl. Acad. Sci. USA* 94:1041–1046.
- Clark J.M. 1988. Novel non-templated nucleotide addition reactions catalyzed by procaryotic and eucaryotic DNA polymerases. *Nucleic Acids Res.* 16:9677–9686.
- Comey, C.T., Koons, B.W., Presley, K.W., Smerick, J.B., Sobieralski, C.A., Stanley, D.M., and Baechtel, F.S. 1994. DNA extraction strategies for amplified fragment length polymorphism analysis. *J. Forensic Sci.* 39:1254–1269.
- Cone, R.W., and Fairfax, M.R. 1993. Protocol for ultraviolet irradiation of surfaces to reduce PCR contamination. *PCR Methods Appl.* 3:S15–S17.
- D2S1338. Cooperative Human Linkage Center (CHLC) accession number 41445. GenBank accession number G08202.
- D16S539. Cooperative Human Linkage Center (CHLC) accession number 715. GenBank accession number G07925.
- D19S433 Cooperative Human Linkage Center (CHLC) accession number 135. GenBank accession number G08036.
- DeFranchis, R., Cross, N.C.P., Foulkes, N.S., and Cox, T.M. 1988. A potent inhibitor of Taq DNA polymerase copurifies with human genomic DNA. *Nucleic Acids Res.* 16:10355.
- DNA Advisory Board, Federal Bureau of Investigation, U.S. Department of Justice. 1998. Quality assurance standards for forensic DNA testing laboratories.
- DNA Recommendations. 1994. Report concerning further recommendations of the DNA Commission of the ISFH regarding PCR-based polymorphisms in STR (short tandem repeat) systems. *Intl. J. Legal Med.* 107:159–160.
- Edwards, A., Civitello, A., Hammond, H., and Caskey, C. 1991. DNA typing and genetic mapping with trimeric and tetrameric tandem repeats. *Am. J. Hum. Genet.* 49:746–756.
- Edwards, A., Hammond, H.A., Lin, J., Caskey, C.T., and Chakraborty, R. 1992. Genetic variation at five trimeric and tetrameric tandem repeat loci in four human population groups. *Genomics* 12:241–253.

- Frégeau, C.J., and Fourney, R.M. 1993. DNA typing with fluorescently tagged short tandem repeats: a sensitive and accurate approach to human identification. *Biotechniques* 15:100–119.
- Gill, P., Urquart, A., Millican, E., Oldroyd, N., Watson, S., Sparkes, R., and Kimpton, C.P. 1996. A new method of STR interpretation using inferential logic—development of a criminal intelligence database. *Int. J. Leg. Med.* 109:14–22.
- 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).
- Guo S.W., and Thompson, E.A. 1992. Performing the exact test of Hardy-Weinberg proportion for multiple alleles. *Biometrics* 48:361–372.
- Hammond, H., Jin, L., Zhong, Y., Caskey, C., and Chakraborty, R. 1994. Evaluation of 13 short tandem repeat loci for use in personal identification applications. *Am J. Hum. Genet.* 55:175–189.
- Hartl, D.L., and Clark, A.G. 1989. *Principles of population genetics*, 2nd edition. Sunderland, MA: Sinauer Associates, Inc.
- Karlin, S., Cameron, E.C., and Williams, P.T. 1981. Sibling and parent offspring correlation estimation with variable family size. *Proc. Natl. Acad. Sci. (USA)* 78:2664–2668.
- Kimpton, C., Walton, A., and Gill, P. 1992. A further tetranucleotide repeat polymorphism in the vWF gene. *Hum. Mol. Genet.* 1:287.
- Kimpton, C.P., Gill, P., Walton, A., Urquhart, A., Millican, E.S., and Adams, M. 1993. Automated DNA profiling employing multiplex amplification of short tandem repeat loci. *PCR Methods Appl.* 3:13–22.
- Kimpton, C.P., Oldroyd, N.J., Watson, S.K., Frasier, R.R.E., Johnson, P.E., Millican, E.S., Urquhart, A., Sparkes, R.L., and Gill, P. 1996. Validation of highly discriminating multiplex short tandem repeat amplification systems for human identification. *Electrophoresis* 17: 1283-1293.
- Kwok, S., and Higuchi, R. 1989. Avoiding false positives with PCR. *Nature* 339:237–238.
- Lazaruk, K., Walsh, P.S., Oaks, F., Gilbert, D., Rosenblum, B.B., Menchen, S., Scheibler, D., Wenz, H.M., Holt, C., Wallin, J. 1998. Genotyping of forensic short tandem repeat (STR) systems based on sizing precision in a capillary electrophoresis instrument. *Electrophoresis* 19:86–93.
- Lareu, M.-V., Barral, S., Salas, A., Rodriguez, M., Pestoni, C., and Carracedo, A. 1998. Further exploration of new STRs of interest for forensic genetic analysis. In: Olaisen, B., Brinkmann, B., and Lincoln, P.J., eds. *Progress in Forensic Genetics 7: Proceedings of the 17th International ISFH Congress, Oslo 2-6 September 1997*. Elsevier, Amsterdam, pp. 192-200.
- Li, H. Schmidt, L., Wei, M-H., Hustad, T. Leman, M.I., Zbar, B., and Tory, K. 1993. Three tetranucleotide polymorphisms for loci:D3S1352; D3S1358; D3S1359. *Hum. Mol. Genet.* 2:1327.
- Luna, L.G., ed. *Manual of Histologic Staining Methods of the Armed Forces Institute of Pathology*, McGraw-Hill Book Co., New York (1968).

- Magnuson, V.L., Ally, D.S., Nylund, S.J., Karanjawala, Z.E., Rayman, J.B., Knapp, J.I., Lowe, A.L., Ghosh, S., and Collins, F.S. 1996. Substrate nucleotide-determined non-templated addition of adenine by Taq DNA polymerase: implications for PCR-based genotyping and cloning. *Biotechniques* 21:700–709.
- Mancuso, D.J., Tuley, E.A., Westfield, L.A., Worrall, N.K., Shelton-Inloes, B.B., Sorace, J.M., Alevy, Y.G., and Sadler, J.E. 1989. Structure of the gene for human von Willebrand factor. *J. Biol. Chem.* 264:19514–19527.
- Meyer, E., Wiegand, P., Rand, S., Kuhlmann, D., Brack, M., and Brinkmann, B. 1995. Microsatellite polymorphisms reveal phylogenetic relationships in primates. *J. Mol. Evol.* 41(1):10–4.
- Mills, K.A., Even, D., and Murrau, J.C. 1992. Tetranucleotide repeat polymorphism at the human alpha fibrinogen locus (FGA). *Hum. Mol. Genet.* 1:779.
- Möller, A., Meyer, E., and Brinkmann, B. 1994. Different types of structural variation in STRs: HumFES/FPS, HumVWA, and HumD21S11. *Intl. J. Legal Med.* 106:319–323.
- Möller, A., and Brinkmann, B. 1995. PCR-VNTRs (PCR—Variable Number of Tandem Repeats) in forensic science. *Cell. Molec. Biol.* 41:715–724.
- Momhinweg, E., Luckenbach, C., Fimmers, R., and Ritter, H. 1998. D3S1358: sequence analysis and gene frequency in a German population. *Forensic Sci. Int.* 95:173–178.
- Nakahori, Y., Takenaka, O., and Nakagome, Y. 1991. A human X-Y homologous region encodes amelogenin. *Genomics* 9:264–269.
- National Research Council. 1996. *The evaluation of forensic DNA evidence*. National Academy Press, Washington, D.C.
- Nei, M., and Roychoudhury, A.K. 1974. Sampling variances of heterozygosity and genetic distance. *Genetics* 76:379–390.
- Nei, M. 1978. Estimation of average heterozygosity and genetic distance from a small number of individuals. *Genetics* 89:583–590.
- Oldroyd, N.J., Urquhart, A.J., Kimpton, C.P., Millican, E.S., Watson, S.K., Downes, T., and Gill, P.D. 1995. A highly discriminating octoplex short tandem repeat polymerase chain reaction system suitable for human individual identification. *Electrophoresis* 16:334–337.
- Oppitz, E. 1969. *Arkiv Fur Kriminologie* 1969:144–145.
- Perkin-Elmer Corporation. 1993. *AmpliType™ User's Guide*, Version 2, Foster City, CA.
- Prince, A.M., and Andrus, L. 1992. PCR—How to kill unwanted DNA. *Biotechniques* 12:358.
- Puers, C., Hammond, H., Jin, L., Caskey, C., and Schumm, J. 1993. Identification of repeat sequence heterogeneity at the polymorphic short tandem repeat locus HUMTH01 [AATG]<sub>n</sub> and reassignment of alleles in population analysis using a locus-specific allelic ladder. *Am. J. Hum. Genet.* 53:953–958.
- Revised Validation Guidelines-Scientific Working Group on DNA Analysis Methods (SWGDM). *Forensic Science Communications* (July 2004) Volume 6 (3). Available at [www.fbi.gov/hq/lab/fsc/current/standards/2004\\_03\\_standards02.htm](http://www.fbi.gov/hq/lab/fsc/current/standards/2004_03_standards02.htm)
- Sambrook, J., Fritsch, E.F., and Maniatis, T., eds. 1989. *Molecular Cloning: A Laboratory Manual, 2nd Edition, Volume 2*. Cold Spring Harbor Laboratory Press, New York, pp. E10–E14.

- Sensabaugh, G.F. 1982. Biochemical markers of individuality. In: Saferstein, R., ed. *Forensic Science Handbook*. Prentice-Hall, Inc., New York, pp. 338–415.
- Sensabaugh G.F., von Beroldingen C. The polymerase chain reaction: application to the analysis of biological evidence. In: Farley, MA, Harrington, JJ, editors. *Forensic DNA Technology*. Michigan: Lewis, 1991:63-82.
- Sharma, V., and Litt, M. 1992. Tetranucleotide repeat polymorphism at the D21S11 locus. *Hum Mol. Genet.* 1:67.
- Singer-Sam, J., and Tanguay, R. 1989. Use of Chelex to Improve the PCR Signal From a Small Number of Cells. *Amplifications* 3.
- Smith, R.N. 1995. Accurate size comparison of short tandem repeat alleles amplified by PCR. *Biotechniques* 18:122–128.
- Sparkes, R., Kimpton, C., Watson, S., Oldroyd, N., Clayton, T., Barnett, L., Arnold, J., Thompson, C., Hale, R., Chapman, J., Urquhart, A., and Gill, P. 1996a. The validation of a 7-locus multiplex STR test for use in forensic casework. (I). Mixtures, ageing, degradation and species studies. *Int. J. Legal Med.* 109:186–194.
- Sparkes, R., Kimpton, C., Gilbard, S., Carne, P., Andersen, J., Oldroyd, N., Thomas, D., Urquhart, A., and Gill, P. 1996b. The validation of a 7-locus multiplex STR test for use in forensic casework. (II), Artifacts, casework studies and success rates. *Int. J. Legal Med.* 109:195–204.
- Straub, R.E., Speer, M.C., Luo, Y., Rojas, K., Overhauser, J., Ott, J., and Gilliam, T.C. 1993. A microsatellite genetic linkage map of human chromosome 18. *Genomics* 15:48–56.
- Sullivan, K.M., Mannucci, A., Kimpton, C.P., and Gill, P. 1993. A rapid and quantitative DNA sex test: fluorescence-based PCR analysis of X-Y homologous gene amelogenin. *Biotechniques* 15:636–641.
- Szibor, R., Lautsch, S., Plate, I., Bender, K., and Krause, D. 1998. Population genetic data of the STR HumD3S1358 in two regions of Germany. *Int. J. Legal Med.* 111:160–161.
- Technical Working Group on DNA Analysis Methods. 1995. Guidelines for a quality assurance program for DNA analysis. *Crime Lab. Digest* 22: 21–43.
- Urquhart, A., Oldroyd, N.J., Kimpton, C.P., and Gill, P. 1995. Highly discriminating heptaplex short tandem repeat PCR system for forensic identification. *Biotechniques* 18:116–121.
- U.S. Department of Health and Human Services. 1993. *Biosafety in Microbiological and Biomedical Laboratories*, 3rd edition, U.S. Government Printing Office.
- U.S. Department of Health and Human Services. OSHA Bloodborne Pathogen Standard 29 CFR, part 1910.1030.
- Wallin, J.M., Buoncristiani, M.R., Lazaruk, K.D., Fildes, N., Holt, C.L., Walsh, P.S. 1998. TWGDAM validation of the AmpFISTR blue PCR amplification kit for forensic casework analysis. *J. Forensic Sci.* 43:854–870.
- Walsh, P.S., Fildes, N.J., and Reynolds, R. 1996. Sequence analysis and characterization of stutter products at the tetranucleotide repeat locus vWA. *Nucleic Acids Res.* 24:2807–2812.

Watson, S., Kelsey, Z., Webb, R., Evans, J., and Gill, P. 1998. The development of a third generation STR multiplex system (TGM). In: Olaisen, B., Brinkmann, B., and Lincoln, P.J., eds. *Progress in Forensic Genetics 7: Proceedings of the 17th International ISFH Congress, Oslo 2-6 September 1997*. Elsevier, Amsterdam, pp. 192–194.

Weber, J. and Wong, C. 1993. Mutation of human short tandem repeats. *Hum. Mol. Genet.* 2:1123–1128.

Weir, B.S. 1996. *Genetic data analysis II*. Sunderland, MA: Sinauer Associates, Inc.

Weir, B.S. 1992. Independence of VNTR alleles defined by fixed bins. *Genetics* 130: 873-887.

Ziegle, J.S., Su, Y., Corcoran, K.P., Nie, L., Mayrand, P.E., Hoff, L.B., McBride, L.J., Kronick, M.N., and Diehl, S.R. 1992. Application of automated DNA sizing technology for genotyping microsatellite loci. *Genomics* 14:1026–1031.

# Documentation and Support

## Related documentation

Document title	Pub. no.
<i>AmpFtSTR™ SGM Plus™ PCR Amplification Kit: Human Identification: Application Note</i>	040302
<i>3100/3100-Avant Data Collection v2.0 User Guide</i>	4347102
<i>3100/3100-Avant Genetic Analyzers Using Data Collection Software v2.0 User Bulletin</i>	4350218
<i>3100 Genetic Analyzer User Manual (Data Collection v1.1)</i>	4315834
<i>3100/3100-Avant Genetic Analyzers Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin</i>	4332345
<i>Applied Biosystems 3130/3100xl Genetic Analyzers Using Data Collection Software v3.0 User Bulletin</i>	4363787
<i>Applied Biosystems 3130/3130xl Genetic Analyzers Getting Started Guide</i>	4352715
<i>Applied Biosystems 3130/3130xl Genetic Analyzers Maintenance, Troubleshooting, and Reference Guide</i>	4352716
<i>Applied Biosystems 3130/3130xl Genetic Analyzers Quick Reference Card</i>	4362825
<i>Applied Biosystems 3130/3130xl Genetic Analyzers AB Navigator Software Administrator Guide</i>	4359472
<i>Applied Biosystems 3130/3100xl DNA Analyzers User Guide</i>	4331468
<i>Applied Biosystems 3500/3500xL Genetic Analyzer Quick Reference Card</i>	4401662
<i>Applied Biosystems 3500/3500xL Genetic Analyzer User Guide, Data Collection v1.0</i>	4401661
<i>Applied Biosystems 3500/3500xL Genetic Analyzer User Bulletin: Solutions to issues related to software, data, hardware, and consumables</i>	4445098
<b>Note:</b> Additional user bulletins may be available at <a href="http://www.lifetechnologies.com">www.lifetechnologies.com</a>	
<i>Applied Biosystems 3730/3730xl Genetic Analyzer Getting Started Guide</i>	4359476
<i>GeneAmp™ PCR System 9700 Base Module User's Manual</i>	N805-0200
<i>Veriti™ 96-Well Thermal Cycler AmpFtSTR™ Kit Validation User Bulletin</i>	4440754
<i>Quantifiler™ Kits: Quantifiler™ Human DNA Quantification Kit and Quantifiler™ Y Human Male DNA Quantification Kit User's Manual</i>	4344790
<i>PrepFiler™ Forensic DNA Extraction Kit User Guide</i>	4390932
<i>GeneMapper™ ID Software Version 3.1 Human Identification Analysis User Guide</i>	4338775
<i>GeneMapper™ ID Software Versions 3.1 and 3.2 Human Identification Analysis Tutorial</i>	4335523
<i>Installation Procedures and New Features for GeneMapper™ ID Software v3.2 User Bulletin</i>	4352543
<i>GeneMapper™ ID-X Software Version 1.0 Getting Started Guide</i>	4375574
<i>GeneMapper™ ID-X Software Version 1.0 Quick Reference Guide</i>	4375670
<i>GeneMapper™ ID-X Software Version 1.0 Reference Guide</i>	4375671
<i>GeneMapper™ ID-X Software Version 1.1 (Mixture Analysis) Getting Started Guide</i>	4396773

Document title	Pub. no.
<i>GeneMapper™ ID-X Software Version 1.1 (Mixture Analysis) Quick Reference Guide</i>	4402094
<i>GeneMapper™ ID-X Software Version 1.2 Reference Guide</i>	4426481
<i>GeneMapper™ ID-X Software Version 1.2 Quick Reference Guide</i>	4426482

Portable document format (PDF) versions of this guide and the documents listed above are available at [www.lifetechnologies.com](http://www.lifetechnologies.com).

**Note:** To open the user documentation available from the Applied Biosystems web site, use the Adobe ▶ Acrobat ▶ Reader ▶ software available from [www.adobe.com](http://www.adobe.com).

## Obtain SDSs

Safety Data Sheets (SDSs) are available from [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

**Note:** For the SDSs of chemicals not distributed by Life Technologies, contact the chemical manufacturer.

## Obtain support

For HID support:

- **In North America** – Send an email to [HIDTechSupport@lifetech.com](mailto:HIDTechSupport@lifetech.com), or call 888-821-4443 option 1.
- **Outside North America** – Contact your local support office.

For the latest services and support information for all locations, go to:

[www.lifetechnologies.com](http://www.lifetechnologies.com)

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant this product as set forth in the *Applied Biosystems™ 3500/3500xL Genetic Analyzer – Instrument Limited Warranty*. If you have any questions, please contact Life Technologies at [thermofisher.com/support](http://thermofisher.com/support).

## A

A nucleotide, addition by AmpliTaq Gold to 3' end of amplicon 77

agarose gel, using to examine DNA 89

allele frequencies in the population databases 93–99

allelic dropout 108

allelic ladder

- about 18
- accuracy 68
- precision 70
- profile 12
- requirements for accurate genotyping 25
- using to determine genotypes 68
- volume per reaction 28, 30, 32

amplification

- differential amplification of loci 89
- loci 11
- using bloodstained FTA cards 23

AmpliTaq Gold DNA Polymerase, catalyzing the addition of a 3' A nucleotide 77

## B

bins

- check version 49
- import 35, 50

biohazard safety 123

buffer, new 102

## C

Caucasian 99

chemical safety 122

contents of kit 18

control DNA 007 14, 18

## D

Data Collection Software, overview 16

developmental validation 66

differential amplification of loci 89

DNA

control, about 18

effect of DNA quantity on results 85

how degraded DNA affects which loci amplify 88, 91

mixed samples 77, 78, 79, 81

negative-control reaction 21

positive-control reaction 21

quantification 19

quantification methods 20

sample preparation 21

test sample 21

using agarose gel analysis to examine the DNA 89

DNA from more than one individual 81

- detecting 78
- detection limit 79, 81

documentation, related 131

## E

electropherogram

- causes for extra peaks 74, 77, 78, 79, 80, 81

electrophoresis

- Data Collection Software 27, 29, 31
- preparing samples on the 310 instrument 31
- preparing samples on the 3100/3100-Avant or 3130/3130xl instrument 28
- preparing samples on the 3500/3500xL instrument 29
- reagents and parts 27, 29, 31
- references 27, 29, 31
- run module 27, 29, 31
- set up 27, 29, 31

emission spectra 17

enzyme, new 102

equipment, not included with kit 117

evidence, exclusion of suspects 92

## F

fluorescent dyes 16

FTA cards

amplification 23  
bloodstained 23

## G

gels 89

GeneMapper ID Software

analyze project 45  
create analysis method 38  
create size standard 43  
examine and edit project 46  
import panels and bins 35  
overview 33  
set up 34

GeneMapper ID-X Software

analyze project 62  
check version of panels, bins, and stutter 49  
create analysis method 55  
create size standard 60  
examine and edit project 63  
import panels, bins, and stutter 50  
overview 48  
set up 49

GeneScan size standard

about 18  
dye label 16  
volume per reaction 28, 29, 31

genetics 92

allele frequencies 93, 99  
populations and samples used in studies 93

genotype

calculating precision data using the allelic ladders 70  
determining 68, 70  
exclusion of suspects 92  
resolving in mixed samples 79

## H

hematin, effect on DNA samples 88

Hi-Di formamide, volume per reaction 28, 29, 31

## I

instrumentation

310 genetic analyzer 16, 31  
3100/3100-Avant genetic analyzer 16, 27  
3130/3130xl genetic analyzer 16, 27  
3500/3500xL genetic analyzer 16, 27, 29  
software compatibility 16

## K

kit

allelic ladder 18  
amplification 11  
contents 18  
control DNA 18  
description 11  
DNA polymerase 18, 21  
fluorescent dyes 16  
loci amplification 11  
PCR reaction mix 18  
primers 11, 18, 20  
reagents 18  
supported instruments 11  
thermal cyclers for use with 116

## L

limited product warranty 133

LIZ size standard

about 18  
volume per reaction 28, 29, 31

loci

allele frequencies in the population  
databases 93–99  
amplified by the kit 11  
chromosomal location 12  
degraded DNA 88, 91  
differential amplification 89  
dye label 12  
effect of DNA quantity on results 85  
effect of inhibitors 88  
lack of amplification 86, 88  
population data 92  
population data, allele frequencies 93  
population data, samples used in studies 93

locus. *See* loci

low-TE buffer 19

## M

master mix, volume per reaction 21

materials and equipment

included in kit 18  
not included with kit 117

multicomponent analysis 16, 17

mutation studies 99

**N**

negative control, sample preparation 21

**O**

operating systems 16, 27, 29, 31

ordering information 117

**P**

panel

check version 49

import 35, 50

part numbers 117

PCR

amplification of tetranucleotide STR loci (stutter peak) 74

cycle number, selecting 22

inhibitor 88

lack of amplification 88

performing 22

setup 115

thermal cycling conditions, programming 22

PCR work areas 115, 117

population genetics 92

allele frequencies 93, 99

populations and samples used in the studies 93

*See Also* allele

population studies 83

positive control, sample preparation 21

precision results 70

primers in the kit 11

primers, volume per reaction 21

probability of identity

definition 100

values 100

product description 11

purpose of kit 11

**Q**

quantification, DNA 19

**R**

reaction mix, for PCR 21

reactions, preparing for PCR 21

reagents

not included with kit 117

user supplied 19

run module, electrophoresis 27, 29, 31

**S**

safety

biohazard 122

chemical 122

Safety Data Sheets (SDSs), obtaining 132

sample preparation 21

DNA negative control 21

DNA positive control 21

standards 18

samples, DNA from more than one individual 77, 81

detecting 78

detection limit 79, 81

resolving genotypes in mixed samples 79

sexual assault

analysis of sexual assault DNA mixtures 82

software, instrument compatibility 16

stutter

check version 49

import 50

stutter peak 74

stutter peak or product 74

support, obtaining 132

**T**

technical support 132

thermal cyclers

for use with kit 116

programming conditions 22

training, information on 132

troubleshooting 113

**U**

user-supplied reagents 19

**V**

validation

developmental 66

importance of 66

mutation rate 99

performance after buffer and enzyme replacement 102

probability of identity 100

probability of paternity exclusion 100

reproducibility [67](#)  
sexual assault DNA mixtures [82](#)

## **W**

warranty [133](#)  
work area  
    amplified DNA [116](#)  
    PCR setup [115](#)  
    setup and lab design [115](#)  
workflow overview [15](#)

