

AmpF ℓ STR[™] NGM[™] PCR Amplification Kit

USER GUIDE

for use with:

200 reaction kit (Part no. 4415020)

1,000 reaction kit (Part no. 4415021)

Catalog Number 4415020, 4415021

Publication Number 4425511

Revision H



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. 4425511

Revision	Date	Description
H	24 August 2018	Updated branding and trademarks, no technical changes.
G	February 2015	Add information for the ProFlex™ PCR System.
F	March 2012	Change copyright information. Added information for GeneAmp™ 9700 and Veriti™ Thermal Cycler emulsion mode in PCR chapter.
E	July 2011	Information on addition of SNP-Specific Primers at D2S441, D22S1045, Amelogenin.
D	July 2011	Patent upgrade.
C	September 2009	Update copyright information.
B	November 2009	Add Chapter 5 Experiments and Results
A	September 2009	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. FTA is a trademark of Whatman Limited. Windows and Windows Vista are trademarks of Microsoft Corporation. Adobe, Acrobat and Reader are trademarks of Adobe Systems Incorporated.

©2018 Thermo Fisher Scientific Inc. All rights reserved.

Contents

	About This Guide	9
	Purpose	9
Chapter 1	Overview	11
	Product overview	11
	Purpose	11
	Product description	11
	About the primers	12
	Loci amplified by the kit	12
	Allelic ladder profile	13
	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	16
	Materials and equipment	17
	Kit contents and storage	17
	Standards for samples	18
Chapter 2	PCR Amplification	19
	Required user-supplied materials and reagents	19
	DNA quantification	19
	Importance of quantification	19
	Methods of quantifying DNA	20
	Prepare the amplification kit reactions	21
	Perform PCR	22
	Amplification using bloodstained FTA™ cards	23
Chapter 3	Electrophoresis	25
	Allelic ladder requirements	25

Section 3.1 3100/3100-*Avant* and 3130/3130*xl* instruments 27
 Set up the 3100/3100-*Avant* and 3130/3130*xl* instruments for electrophoresis 27
 Reagents and parts 27
 Electrophoresis software setup and reference documents 27
 Prepare samples for electrophoresis on the 3100/3100-*Avant* or
 3130/3130*xl* instruments 28

Section 3.2 3500/3500*xL* instruments 29
 Set up the 3500/3500*xL* instrument for electrophoresis 29
 Reagents and parts 29
 Electrophoresis software setup and reference documents 29
 Prepare samples for electrophoresis on the 3500/3500*xL* instrument 29

Section 3.3 310 instruments 31
 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

Chapter 4 Data Analysis 33

Section 4.1 GeneMapper™ *ID* Software 33
 Overview of GeneMapper™ *ID* Software 33
 Instruments 33
 Before you start 34
 Set up GeneMapper™ *ID* Software for data analysis 34
 File names 34
 Overview 34
 Import panels and bins 35
 Create an analysis method 38
 General tab settings 38
 Allele tab settings 39
 Peak Detector tab settings 40
 Peak Quality tab settings 41
 Quality Flags tab settings 42
 Create a size standard 42
 Analyze and edit sample files with GeneMapper™ *ID* Software 43
 Examine and edit a project 44
 For more information 44

Section 4.2 GeneMapper™ *ID-X* Software 45
 Overview of GeneMapper™ *ID-X* Software 45
 Instruments 45
 Before you start 45

Set up GeneMapper™ <i>ID-X</i> Software for data analysis	46
File names	46
Overview	46
Import panels, bins, and marker stutter	46
Create an analysis method	50
General tab settings	51
Allele tab settings	52
Peak Detector tab settings	53
Peak Quality tab settings	54
SQ & GQ tab settings	55
Create a size standard	55
Analyze and edit sample files with GeneMapper™ <i>ID-X</i> Software	57
Examine and edit a project	58
For more information	58
Chapter 5	
Experiments and Results	59
Overview	59
Importance of validation	59
Experiment conditions	59
Developmental validation	60
SWGDM guideline 1.2.1	60
SWGDM guideline 2.10.1	60
PCR components	60
Thermal cycler parameters	61
PCR cycle number	62
Accuracy, precision, and reproducibility	63
SWGDM guideline 2.9	63
Accuracy	63
Precision and size windows	64
Extra peaks in the electropherogram	72
Causes of extra peaks	72
Stutter products	72
Addition of 3' A nucleotide	76
Artifacts	77
Characterization of loci	79
SWGDM guideline 2.1	79
Nature of the polymorphisms	79
Inheritance	80
Mapping	80
Genetic linkage	80
Species specificity	81
SWGDM Guideline 2.2	81

Sensitivity	82
SWGDM guideline 2.3	82
Importance of quantification	82
Effect of DNA quantity on results	83
Stability	84
SWGDM guideline 2.4	84
Degraded DNA	84
Effect of inhibitors — hematin	85
Effect of inhibitors — humic acid	87
Mixture studies	88
SWGDM guideline 2.8	88
Mixture studies	88
Resolution of genotypes in mixed samples	89
Limit of detection of the minor component	89
Population data	91
SWGDM guideline 2.7	91
Overview	91
Population samples used in these studies	91
New Primers added to the NGM™ Kit	91
Comparison of primer sequences with other AmpF \mathcal{L} STR™ Kits	91
Inclusion of three SNP-specific primers to address mutations at the Amelogenin, D2S441, and D22S1045 loci	92
Mutation rate	93
Estimating germ-line mutations	93
Additional mutation studies	94
Probability of identity	94
Probability of paternity exclusion	104
Appendix A Troubleshooting	105
Appendix B Ordering Information	107
Materials and equipment not included	107
Appendix C PCR Work Areas	111
Work area setup and lab design	111
PCR setup work area	111
Amplified DNA work area	112
Appendix D Safety	113
Chemical safety	114

Biological hazard safety 114

Documentation and Support 117

 Related documentation 117

 Obtain support 118

 Limited Product Warranty 118

Bibliography 119

Index 123

About This Guide

Purpose

The *Applied Biosystems™ AmpFℓSTR™ NGM™ PCR Amplification Kit User Guide* provides information about our instruments, chemistries, and software associated with the AmpFℓSTR™ NGM™ PCR Amplification Kit.

- Product overview 11
- Workflow overview 15
- Instrument and software overview 16
- Materials and equipment 17

Product overview

Purpose

The AmpF ℓ STR[™] NGM[™] PCR Amplification Kit is a short tandem repeat (STR) multiplex assay that amplifies 14 tetranucleotide repeat loci and one trinucleotide repeat locus, D22S1045. The kit simultaneously coamplifies the 10 loci contained in the AmpF ℓ STR[™] SGM Plus[™] Kit (D3S1358, vWA, D16S539, D2S1338, D8S1179, D21S11, D18S51, D19S433, TH01, and FGA) together with 2 highly polymorphic STR loci (D1S1656 and D12S391), 3 “mini” STR loci (D10S1248, D22S1045 and D2S441), and the gender determination locus Amelogenin. The NGM[™] Kit delivers a 16-locus multiplex with a greater power of discrimination, better sensitivity, and improved robustness than earlier generation AmpF ℓ STR[™] kits. The kit uses modified PCR cycling conditions for enhanced sensitivity, a new buffer formulation to improve performance with inhibited samples, more loci concentrated in the low molecular-weight region of the profile to improve performance on degraded samples, and an improved process for synthesis and purification of the amplification primers to deliver a much cleaner electrophoretic background.

Product description

The NGM[™] Kit contains all the necessary reagents for the amplification of human genomic DNA.

The reagents are designed for use with the following Applied Biosystems[™] instruments:

- ABI PRISM[™] 3100/3100-*Avant* Genetic Analyzer
- Applied Biosystems[™] 3130/3130*xl* Genetic Analyzer
- Applied Biosystems[™] 3500/3500*xL* Genetic Analyzer
- Applied Biosystems[™] 310 Genetic Analyzer
- GeneAmp[™] PCR System 9700 with the Silver 96-Well Block
- GeneAmp[™] PCR System 9700 with the Gold-plated Silver 96-Well Block
- Veriti[™] 96-Well Thermal Cycler
- ProFlex[™] PCR System

About the primers

The NGM™ Kit employs the latest improvements in primer synthesis and purification techniques to minimize the presence of dye-labeled artifacts. These improvements result in a much cleaner electropherogram background that enhances the assay's signal-to-noise ratio and simplifies the interpretation of results.

Loci amplified by the kit

The following table shows the loci amplified, their chromosomal locations, and the corresponding fluorescent marker dyes. The AmpF λ STR™ NGM™ Allelic Ladder is used to genotype the analyzed samples. The alleles contained in the allelic ladder and the genotype of the AmpF λ STR™ Control DNA 007 are also listed in the table.

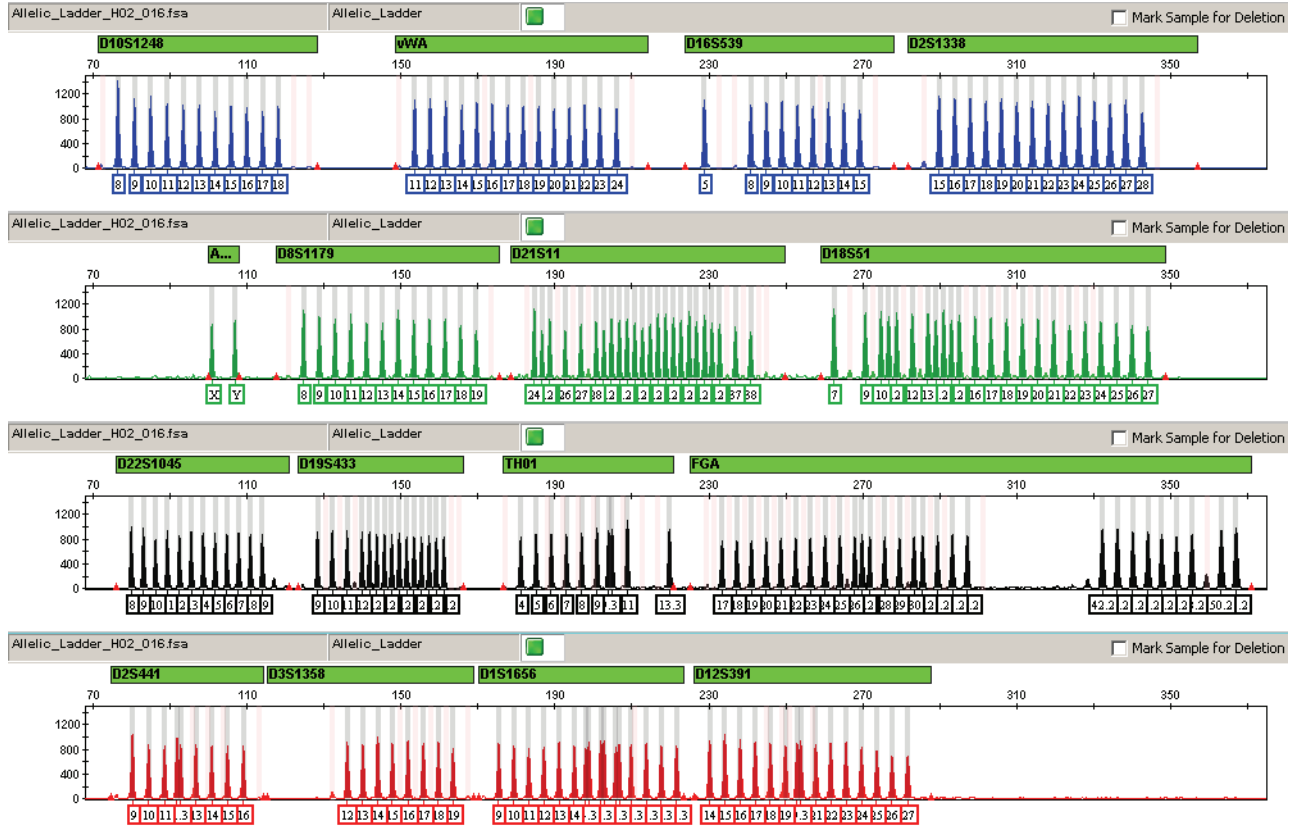
Table 1 NGM™ Kit loci and alleles

Locus designation	Chromosome location	Alleles included in AmpF λ STR™ NGM™ Allelic Ladder	Dye label	Control DNA 007
D10S1248	10q26.3	8, 9, 10, 11, 12, 13, 14, 15, 16, 17, 18	6-FAM™	12, 15
vWA	12p13.31	11,12, 13, 14, 15, 16, 17, 18, 19, 20, 21, 22, 23, 24	6-FAM™	14, 16
D16S539	16q24.1	5, 8, 9, 10, 11, 12,13, 14, 15	6-FAM™	9, 10
D2S1338	2q35	15, 16, 17, 18, 19, 20, 21, 22, 23, 24, 25, 26, 27, 28	6-FAM™	20, 23
Amelogenin	X: p22.1-22.3 Y: p11.2	X, Y	VIC™	X, Y
D8S1179	8q24.13	8, 9 10, 11, 12, 13, 14, 15, 16, 17, 18, 19	VIC™	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	VIC™	28, 31
D18S51	18q21.33	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	VIC™	12, 15
D22S1045	22q12.3	8, 9, 10, 11, 12, 13, 14, 15, 16, 17, 18, 19	NED™	11, 16
D19S433	19q12	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	NED™	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	NED™	24, 26
D2S441	2p14	9, 10, 11, 11.3, 12, 13, 14, 15, 16	PET™	14, 15
D3S1358	3p21.31	12, 13, 14, 15, 16, 17, 18, 19	PET™	15, 16
D1S1656	1q42.2	9, 10, 11, 12, 13, 14, 14.3, 15, 15.3, 16, 16.3, 17, 17.3, 18.3, 19.3, 20.3	PET™	13, 16
D12S391	12p13.2	14, 15, 16, 17, 18, 19, 19.3, 20, 21, 22, 23, 24, 25, 26, 27	PET™	18, 19

Allelic ladder profile

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

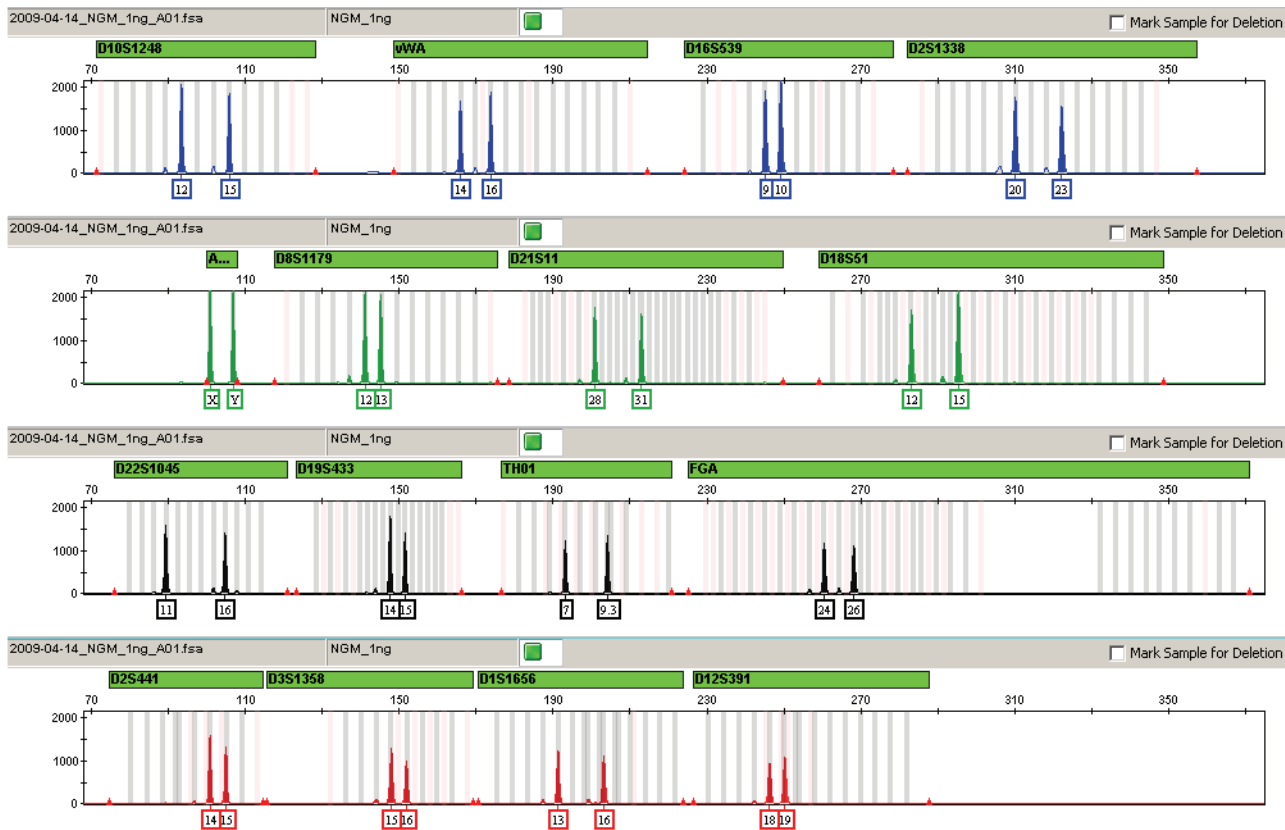
Figure 1 GeneMapper™ /D-X Software plot of the AmpF/STR™ NGM™ Allelic Ladder



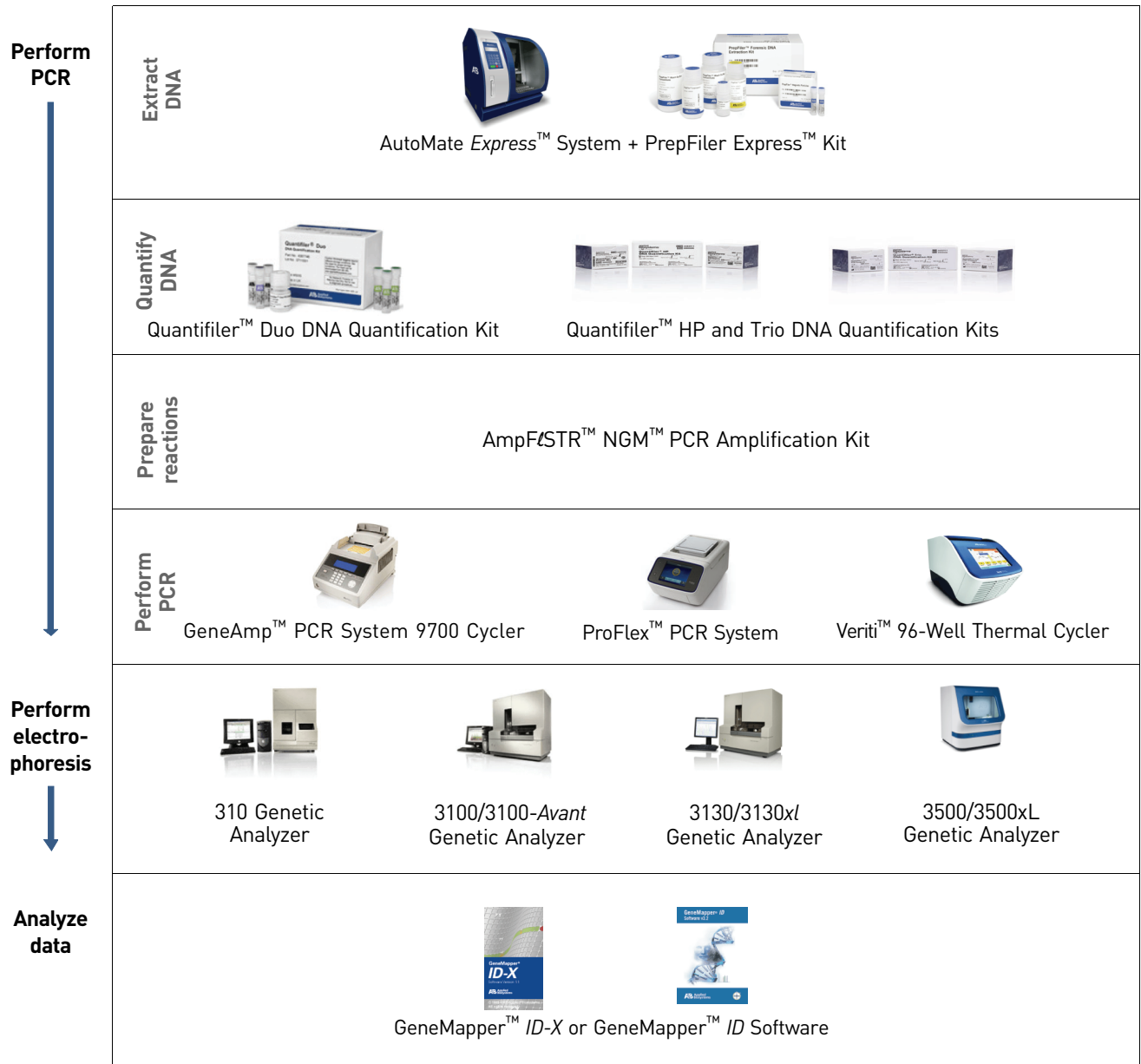
Control DNA 007 profile

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

Figure 2 1 ng of Control DNA 007 amplified with the NGM™ 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 NGM™ 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 files for 31xx instruments and .hid files for 3500 instruments), which is then analyzed by the GeneMapper™ ID or ID-X Software.

Instrument and software compatibility

Table 2 Software specific to each instrument

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

† We conducted validation studies for the NGM™ Kit using this configuration.

About multicomponent analysis

Our 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.

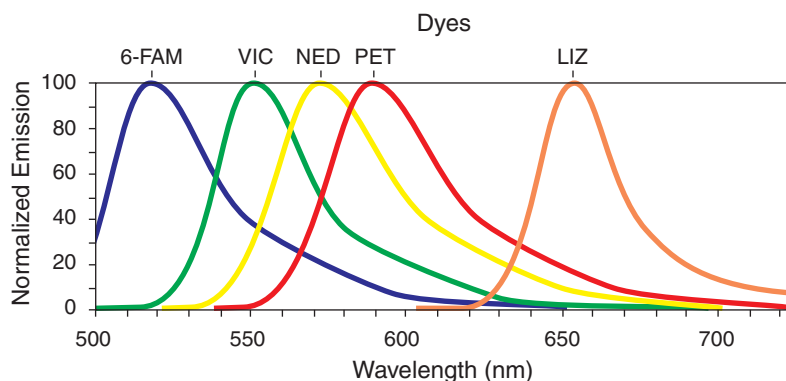
Multicomponent analysis is the process that separates the 5 different fluorescent dye colors into distinct spectral components. The four dyes used in the NGM™ Kit to label samples are 6-FAM™, VIC™, NED™, and PET™ dyes. The fifth dye, LIZ™, is used to label the GeneScan™ 500 LIZ™ Size Standard or GeneScan™ 600 LIZ™ Size Standard v2.0.

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 ABI PRISM™ 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 6-FAM™ dye emits at the shortest wavelength and it is displayed as blue, followed by the VIC™ dye (green), NED™ dye (yellow), PET™ dye (red), and LIZ™ dye (orange).

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 five dyes used in the NGM™ Kit



Materials and equipment

Kit contents and storage

The NGM™ Kit contains materials sufficient to perform 200 (Part no. 4415020) or 1000 (Part no. 4415021) amplifications at a 25 μ L reaction volume.

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

Table 3 Kit contents and storage

Component	Description	200× Volume	1000× Volume	Storage
AmpF λ STR™ NGM™ Primer Set	Contains forward and reverse primers to amplify human DNA targets.	1 tube, 1.0 mL	1 bottle, 5.0 mL	-15 to -25°C on receipt, 2 to 8 °C after initial use
AmpF λ STR™ NGM™ Master Mix	Contains enzyme, salts, dNTPs, carrier protein, and 0.05% sodium azide.	2 tubes, 1.0 mL each	1 bottle, 10.0 mL	-15 to -25°C on receipt, 2 to 8 °C after initial use
AmpF λ STR™ NGM™ Allelic Ladder	Contains amplified alleles. See Table 1 on page 12 for a list of alleles included in the allelic ladder.	1 tube, 50.0 μ L	1 tube, 75.0 μ L	-15 to -25°C on receipt, 2 to 8°C after initial use
AmpF λ STR™ Control DNA 007	Contains 0.10 ng/ μ L human male 007 DNA in 0.02% sodium azide and buffer†. See Table 1 on page 12 for profile.	1 tube, 0.3 mL	1 tube, 0.3 mL	2 to 8°C

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

Standards for samples

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

- **Control DNA 007** – A positive control for evaluating the efficiency of the amplification step and STR genotyping using the AmpF ϕ STR™ NGM™ Allelic Ladder.
- **GeneScan™ 500 LIZ™ Size Standard (GS 500) or GeneScan™ 600 LIZ™ Size Standard v2.0 (GS 600 v2.0)** – Used for obtaining sizing results. These standards, which have been evaluated as internal size standards, yield precise sizing results for NGM™ Kit PCR products. Order the GeneScan™ 500 LIZ™ Size Standard (Part no. 4322682) or the GeneScan™ 600 LIZ™ Size Standard v2.0 (Part no. 4408399) separately.
- **AmpF ϕ STR™ NGM™ Allelic Ladder** – Allelic ladder for accurate characterization of the alleles amplified by the NGM™ Kit. The AmpF ϕ STR™ NGM™ Allelic Ladder contains most of the alleles reported for the 15 autosomal loci. Refer to Table 1 on page 12 for a list of the alleles included in the AmpF ϕ STR™ NGM™ Allelic Ladder.

■ Required user-supplied materials and reagents	19
■ DNA quantification	19
■ Prepare the amplification kit reactions	21
■ Perform PCR	22
■ Amplification using bloodstained FTA™ cards	23

Required user-supplied materials and reagents

In addition to the NGM™ 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. no. 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 based on your 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 NGM™ Kit is 1.0 ng in a maximum input volume of 10 µL amplified for 29 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:
 - Quantification (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

We provide several kits for quantifying DNA in samples. See the references 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> (Part no. 4344790)	<p>Properties: The Quantifiler™ Human and Quantifiler™ Y Human Male Kits are highly specific for human DNA, and they detect total human or male DNA, respectively. The kits detect single-stranded and degraded DNA.</p> <p>How they work: The Quantifiler™ DNA Quantification Kits consist of target-specific and internal control 5' nuclease assays. 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 that consists of an IPC template DNA (a synthetic sequence not found in nature), two primers for amplifying the IPC template DNA, and one TaqMan™ MGB probe labeled with VIC™ dye for detecting the amplified IPC DNA.</p>
Quantifiler™ Duo DNA Quantification Kit (Part no. 4387746) For more information, see <i>Quantifiler™ Duo DNA Quantification Kit User's Manual</i> (Part no. 4391294)	<p>Properties: The Quantifiler™ Duo Kit is highly specific for human DNA and combines the detection of both total human and male DNA in one PCR reaction. The kit detects single-stranded and degraded DNA.</p> <p>How it works: The Quantifiler™ Duo DNA Quantification Kit consists of target-specific and internal control 5' nuclease assays. 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>

Product	Description
<p>Quantifiler™ HP DNA Quantification Kit (Cat. no. 4482911)</p> <p>Quantifiler™ Trio DNA Quantification Kit (Cat. no. 4482910)</p> <p>For more information, see <i>Quantifiler HP and Trio DNA Quantification Kits User Guide</i> (Pub no. 4485354)</p>	<p>Properties:</p> <p>The Quantifiler™ HP Kit is designed to quantify the total amount of amplifiable human DNA in a sample. The Quantifiler™ Trio Kit is designed to simultaneously quantify the total amount of amplifiable human DNA and human male DNA in a sample.</p> <p>How they work:</p> <p>The Quantifiler™ HP and Trio DNA Quantification Kits use multiple-copy target loci for improved detection sensitivity. The human-specific target loci (Small Autosomal, Large Autosomal, and Y-chromosome targets) each consist of multiple copies dispersed on various autosomal chromosomes (Small Autosomal and Large Autosomal).</p> <p>To maximize the consistency of quantification results, genomic targets were selected with conserved primer- and probe-binding sites within individual genomes and also with minimal copy number variability between different individuals and population groups. As a result, the detection sensitivity of the Quantifiler™ HP and Trio assays is improved over Quantifiler™ Duo, Human, and Y Human Male DNA Quantification Kit assays. The primary quantification targets (Small Autosomal and Y) consist of relatively short amplicons (75 to 80 bases) to improve the detection of degraded DNA samples. In addition, the Quantifiler™ HP and Trio Kits each contain a Large Autosomal target with a longer amplicon (>200 bases) to aid in determining if a DNA sample is degraded.</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 ℓ STR™ NGM™ Master Mix	10.0 μ L
AmpF ℓ STR™ NGM™ Primer Set	5.0 μ L

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

2. Prepare reagents. Thaw the AmpF ℓ STR™ NGM™ Master Mix and the AmpF ℓ STR™ NGM™ Primer Set, then vortex the tubes for 3 seconds and centrifuge them briefly before opening.

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

3. Pipet the required volumes of components into an appropriately sized polypropylene tube.
4. Vortex the reaction mix for 3 seconds, then centrifuge briefly.
5. Dispense 15 μ L of 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 10 μL 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 ng of total DNA is in a final volume of 10 μL . Add 10 μL of the diluted sample to the reaction mix.
Positive control	Add 10 μL of 007 control DNA (0.1 ng/ μL) to provide 1.0 ng of total DNA in the positive control reaction.

The final reaction volume (sample or control plus reaction mix) should be 25 μL .

- Seal the MicroAmp™ Optical 96-Well Reaction Plate with MicroAmp™ Clear Adhesive Film or MicroAmp™ Optical Adhesive Film, or cap the tubes.
- Centrifuge the tubes or plate at 3000 rpm for about 20 seconds in a tabletop centrifuge (with plate holders if using 96-well plates) to remove bubbles.
- Amplify the samples in a GeneAmp™ PCR System 9700 with the Silver 96-well block, or a GeneAmp™ PCR System 9700 with the Gold-plated Silver 96-well block, or a Veriti™ 96-well Thermal Cycler, or a ProFlex™ PCR System.

IMPORTANT! The NGM™ 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 the performance of the NGM™ Kit.

Perform PCR

- 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 AmpFtSTR™ Kit Validation* (Part no. 4440754).
 - When using the ProFlex™ PCR System, refer to the *ProFlex™ PCR System Kit Validation User Bulletin* (Pub. no. 100031595) for more information.

Initial incubation step	Cycle (29/30 cycles)		Final extension	Final hold
	Denature	Anneal		
HOLD	CYCLE		HOLD	HOLD
95°C 11 min	94°C 20 sec	59°C 3 min	60°C 10 min	4°C ∞

IMPORTANT! The NGM™ Kit is validated for use at both 29 and 30 cycles. The optimum conditions for the NGM™ Kit are 29 cycles of amplification with a 1 ng input DNA concentration. If using the NGM™ Kit at 30 cycles, reduce the input DNA concentration to 500 pg. Perform internal validation studies to evaluate kit performance at each cycle number intended for operational use.

2. Load the plate or tubes 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.

3. Start the run.
4. 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

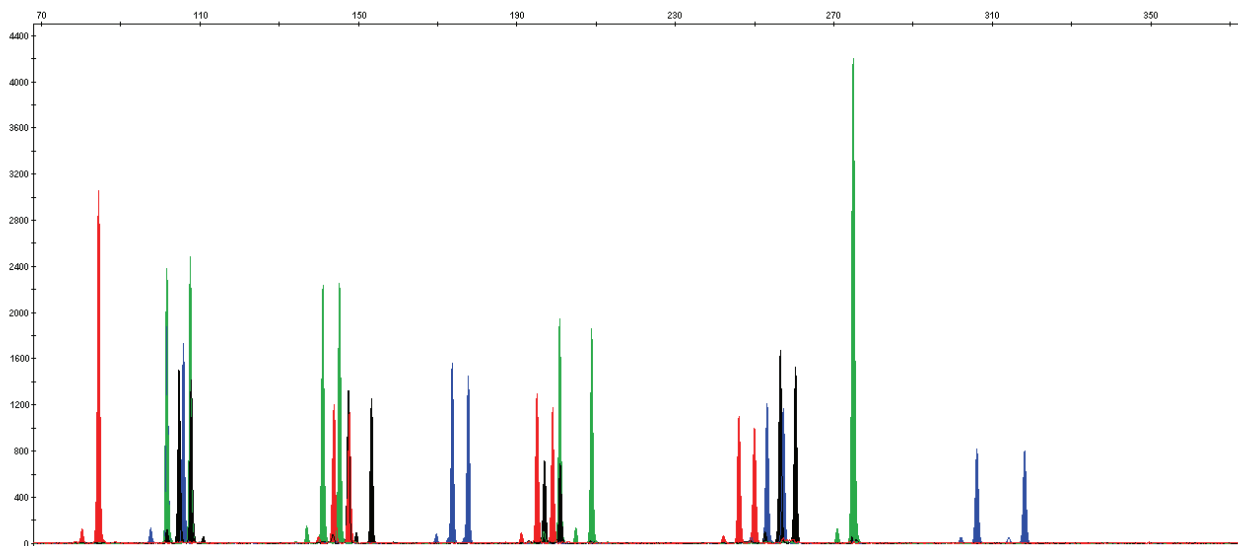
IMPORTANT! Store the amplified products so that they are protected from light.

Amplification using bloodstained FTA™ cards

FTA™ cards can be useful for the collection, storage, and processing of 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 24 cycles, determined by our validation studies. Perform internal validation studies to evaluate kit performance at each cycle number intended for operational use.

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 punch was then amplified directly in the MicroAmp™ tube for 24 cycles.

Figure 4 NGM™ Kit results from a 1.2-mm FTA™ bloodstain disc (24-cycle amplification), analyzed on the Applied Biosystems™ 3130xl Genetic Analyzer



■ Allelic ladder requirements	25
Section 3.1 3100/3100-Avant and 3130/3130xl instruments	27
■ Set up the 3100/3100-Avant and 3130/3130xl instruments for electrophoresis . .	27
■ Prepare samples for electrophoresis on the 3100/3100-Avant or 3130/3130xl instruments	28
Section 3.2 3500/3500xL instruments	29
■ Set up the 3500/3500xL instrument for electrophoresis	29
■ Prepare samples for electrophoresis on the 3500/3500xL instrument.	29
Section 3.3 310 instruments	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 samples	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 and 3130/3130xl instruments for electrophoresis

Reagents and parts Appendix B, “Ordering Information” on page 107 lists the required materials not supplied with the NGM™ 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 NGM™ 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"> HIDFragmentAnalysis36_POP4_1 Injection conditions: <ul style="list-style-type: none"> 3130 = 3 kV/5 sec 3130xl = 3 kV/10 sec Dye Set G5 	<i>Applied Biosystems™ 3130/3130xl Genetic Analyzers Using Data Collection Software v3.0, Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin (Part no. 4363787)</i>
ABI PRISM™ 3100	2.0	Windows™ 2000	<ul style="list-style-type: none"> HIDFragmentAnalysis36_POP4_1 Injection condition: 3kV/10 sec Dye Set G5 	<i>ABI PRISM™ 3100/3100-Avant Genetic Analyzers Using Data Collection Software v2.0, Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin (Part no. 4350218)</i>
	1.1	Windows™ NT	<ul style="list-style-type: none"> GeneScan36vb_DyeSetG5Module Injection condition: 3kV/10 sec GS600v2.0Analysis.gsp 	<i>ABI PRISM™ 3100/3100-Avant Genetic Analyzers Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin (Part no. 4332345)</i>
ABI PRISM™ 3100-Avant	1.0	Windows™ NT	<ul style="list-style-type: none"> GeneScan36Avb_DyeSetG5Module Injection condition: 3 kV/5sec GS600v2.0Analysis.gsp 	<i>ABI PRISM™ 3100/3100-Avant Genetic Analyzers Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin (Part no. 4332345)</i>

† We conducted validation studies for the <Short form of the primary product name> using this configuration.

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

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

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

Reagent	Volume per reaction
GS 500 LIZ™ Size Standard <i>or</i> GS 600 LIZ™ Size Standard v2.0	0.5 µL
Hi-Di™ Formamide	9.5 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. Pipet 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:
 - 10 µL of the formamide:size standard mixture
 - 1 µL of PCR product or Allelic Ladder

Note: For blank wells, add 11 µL of Hi-Di™ Formamide.
5. Seal the reaction plate with appropriate septa, then briefly vortex and centrifuge the plate to ensure that the contents of each well are mixed and 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 on the autosampler.
9. Start the electrophoresis run.

Section 3.2 3500/3500xL instruments

Set up the 3500/3500xL instrument for electrophoresis

Reagents and parts Appendix B, "Ordering Information" on page 107 lists the required materials not supplied with the NGM™ 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 NGM™ 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™ XP	<ul style="list-style-type: none"> HID36_POP4 Injection conditions: 1.2kV/15 sec Dye Set G5 	<i>Applied Biosystems™ 3500/3500xL Genetic Analyzer User Guide</i> (Part no. 4401661) <i>3500 and 3500xL Genetic Analyzers Quick Reference Card</i> (Part no. 4401662)
Applied Biosystems™ 3500xL		Windows Vista™	<ul style="list-style-type: none"> HID36_POP4 Injection conditions: 1.2kV/24 sec Dye Set G5 	

Prepare samples for electrophoresis on the 3500/3500xL instrument

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

1. Calculate the volume of Hi-Di™ Formamide and Size Standard needed to prepare the samples, using the table below.

Reagent	Volume per reaction
GeneScan™ 600 LIZ™ Size Standard v2.0	0.5 µL
Hi-Di™ Formamide	9.5 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. Pipet 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:
 - a. 10 μ L of the formamide: size standard mixture
 - b. 1 μ L of PCR product or allelic ladder
- Note:** For blank wells, add 11 μ L 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 put it onto 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 instruments

Set up the 310 instrument for electrophoresis

Reagents and parts Appendix B, “Ordering Information” on page 107 lists the required materials not supplied with the NGM™ 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 NGM™ 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 [†] or 3.0 [†]	Windows XP or Windows NT™ and Windows 2000	<ul style="list-style-type: none"> GS STR POP4 (1mL) G5 v2.md5 Injection condition: 15 kV/5 sec	ABI PRISM™ 310 Genetic Analyzer User's Manual (Windows) (Part no. 4317588) ABI PRISM™ 310 Protocols for Processing AmpFtSTR™ PCR Amplification Kit Products with Microsoft Windows NT Operating System: User Bulletin (Part no. 4341742)

[†] We conducted concordance studies for the NGM™ Kit using this configuration.

Prepare samples for electrophoresis on the 310 instrument

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

1. Calculate the volume of Hi-Di™ Formamide and Size Standard needed to prepare the samples, using the table below.

Reagent	Volume per reaction
GeneScan™ 500 LIZ™ Size Standard <i>or</i> GeneScan™ 600 LIZ™ Size Standard v2.0	0.75 µL
Hi-Di™ Formamide	24.25 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. Pipet 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:
 - a. 25 μ L of the formamide: size standard mixture
 - b. 1.5 μ L of PCR product or allelic ladder

Note: For blank wells, add 25 μ 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.

Section 4.1 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.....	43
■ Examine and edit a project	44
■ For more information.....	44
Section 4.2 GeneMapper™ ID-X Software	45
■ Overview of GeneMapper™ ID-X Software	45
■ Set up GeneMapper™ ID-X Software for data analysis.....	46
■ Analyze and edit sample files with GeneMapper™ ID-X Software.....	57
■ Examine and edit a project	58
■ For more information.....	58

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 an .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. Perform the appropriate internal validation studies if you want to use multiple ladder samples in an analysis.
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 ± 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, verify the sample result according to your 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.appliedbiosystems.com.

Overview

To analyze sample (.fsa) files using GeneMapper™ ID Software v3.2.1 for the first time:

1. Import panels and bins into the Panel Manager, as explained in “Import panels and bins” on page 35.
2. Create an analysis method, as explained in “Create an analysis method” on page 38.
3. Create a size standard, as explained in “Create a size standard” on page 42.
4. Define custom views of analysis tables.
Refer to the *GeneMapper™ ID Software Versions 3.1 and 3.2 Human Identification Analysis Tutorial* (Part no. 4335523) for more information.
5. Define custom views of plots.
Refer to the *GeneMapper™ ID Software Versions 3.1 and 3.2 Human Identification Analysis Tutorial* (Part no. 4335523) for more information.

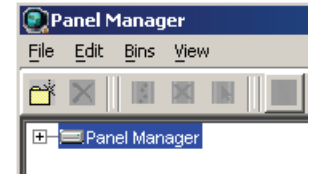
Import panels and bins

To import the NGM™ Kit panel and bin set from www.appliedbiosystems.com into the GeneMapper™ ID Software v3.2.1 database:

1. Download and open the file containing panels and bins:
 - a. From the Support menu of www.appliedbiosystems.com, select **Support ▶ Software Downloads, Patches & Updates ▶ GeneMapper™ ID Software v 3.2 ▶ Updates & Patches**, and download the file **NGM Analysis Files GMID**.
 - b. Unzip the file.
2. Start the GeneMapper™ ID Software, then log in with the appropriate user name and password.

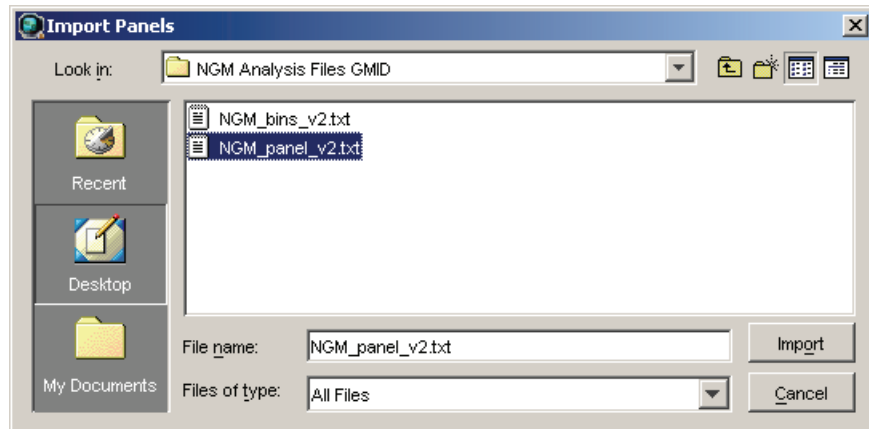
IMPORTANT! For logon instructions, refer to the *GeneMapper™ ID Software Version 3.1 Human Identification Analysis User Guide* (Part no. 4338775).

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.



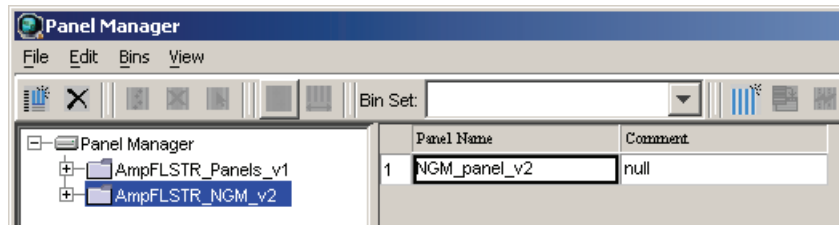
- b. Select **File ▶ Import Panels** to open the Import Panels dialog box.
 - c. Navigate to, then open the **NGM Analysis Files GMID** folder that you unzipped in step 1 above.
5. Select **NGM_panel_v2.txt**, then click **Import**.

Note: Importing this file creates a new folder in the navigation pane of the Panel Manager, AmpFLSTR_NGM_v2. This folder contains the panel and associated markers.



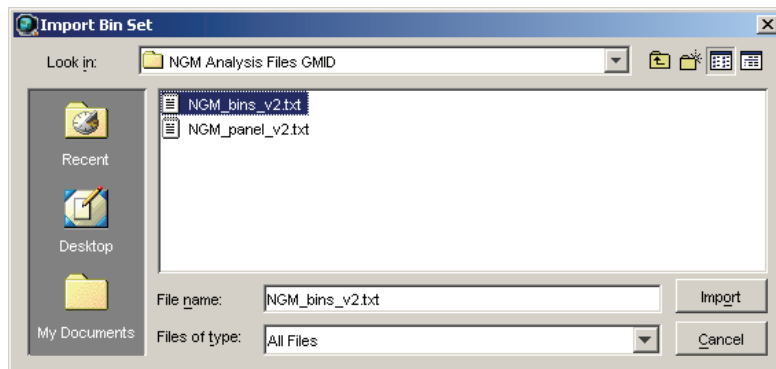
6. Import **NGM_bins_v2.txt**:

- a. Select the **AmpFLSTR_NGM_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 **NGM Analysis Files GMID** folder.
 d. Select **NGM_bins_v2.txt**, then click **Import**.

Note: Importing this file associates the bin set with the panels in the NGM_panel_v2 folder.

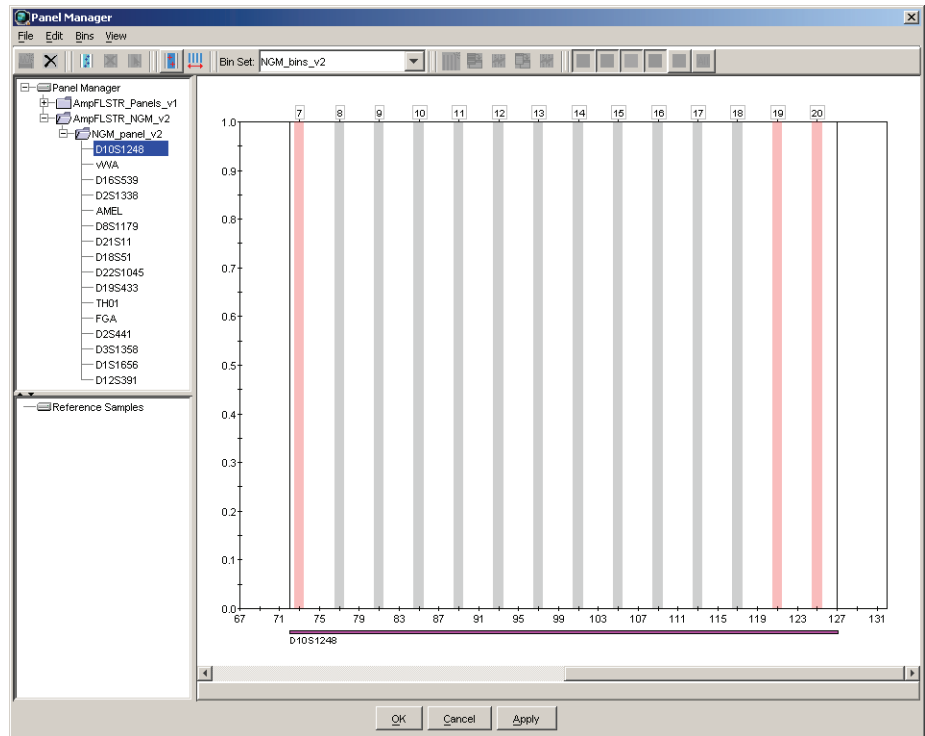


7. View the imported panels in the navigation pane:

- a. Double-click the **AmpFLSTR_NGM_v2** folder to view the **NGM_panel_v2** folder.
 b. Double-click the **NGM_panel_v2** folder to display the panel information in the right pane.

Marker Name	Dye Color	Min Size	Max Size	Control Alleles	Marker I	Marker C	Comments	Ladder Alleles
1 D10S1248	blue	72.0	127.0	"12,15"	4	0.1239	none	8,9,10,11,12,13,14,15,16,17
2 WWA	blue	149.0	214.3	"14,16"	4	0.1183	none	11,12,13,14,15,16,17,18,19,
3 D16S539	blue	223.6	277.6	"9,10"	4	0.1012	none	5,8,9,10,11,12,13,14,15
4 D2S1338	blue	281.6	356.0	"20,23"	4	0.1283	none	15,16,17,18,19,20,21,22,23,
5 AMEL	green	100.0	108.0	"x,y"	9	0.0	none	X,Y
6 D8S1179	green	117.9	174.9	"12,13"	4	0.1031	none	8,9,10,11,12,13,14,15,16,17
7 D21S11	green	178.8	249.8	"28,31"	4	0.1087	none	24,24.2,25,26,27,28,28.2,29
8 D18S51	green	259.5	347.5	"12,15"	4	0.1408	none	7,8,10,10.2,11,12,13,13.2,14
9 D22S1045	yellow	76.0	120.0	"11,16"	3	0.1805	none	8,9,10,11,12,13,14,15,16,17
10 D19S433	yellow	122.3	166.3	"14,15"	4	0.112	none	9,10,11,12,12.2,13,13.2,14,
11 TH01	yellow	176.4	221.1	"7,9,3"	4	0.0427	none	4,5,6,7,8,9,9.3,10,11,13.3
12 FGA	yellow	221.6	372.0	"24,26"	4	0.121	none	17,18,19,20,21,22,23,24,25,
13 D2S441	red	74.5	113.4	"14,15"	4	0.0945	none	9,10,11,11.3,12,13,14,15,16
14 D3S1358	red	114.4	168.4	"15,16"	4	0.1307	none	12,13,14,15,16,17,18,19
15 D1S1656	red	170.0	224.0	"13,16"	4	0.1446	none	9,10,11,12,13,14,14.3,15,15
16 D12S391	red	225.0	287.0	"18,19"	4	0.1527	none	14,15,16,17,18,19,19.3,20,2

8. Select **D10S1248** to display the Bin view for the marker in the right pane.



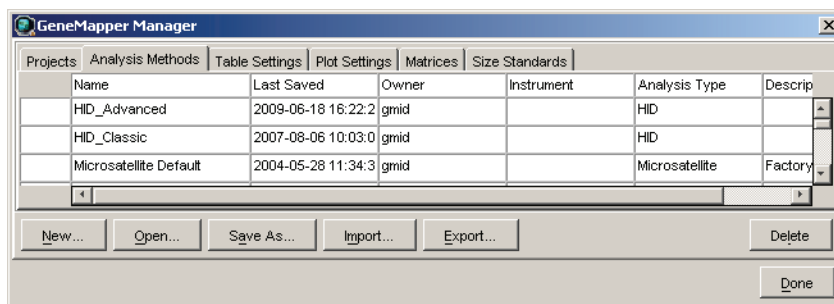
9. Click **Apply**, then **OK** to add the NGM™ Kit panel and bin set to the GeneMapper™ ID Software database.

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

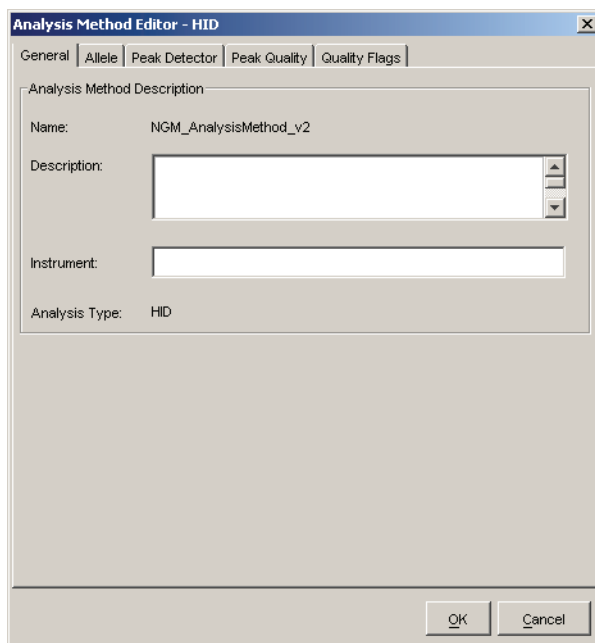
Use the following procedure to create an HID analysis method for the AmpF Φ STR™ NGM™ Kit.

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



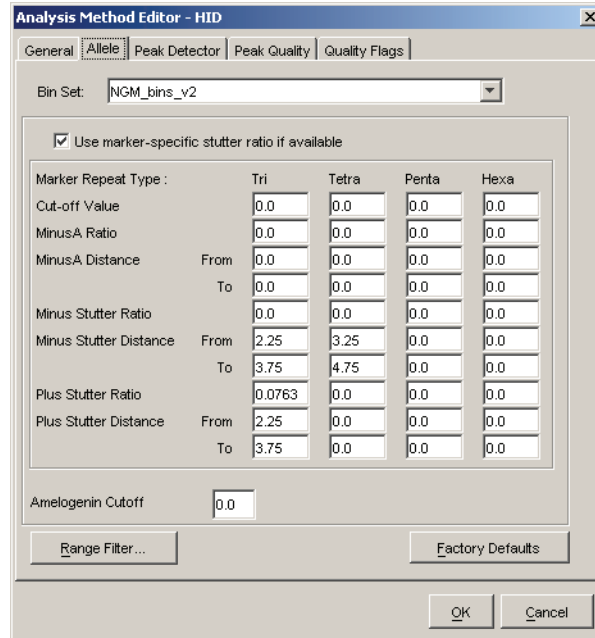
2. Select the **Analysis Methods** tab, then click **New** to open the New Analysis Method dialog box.
3. Select **HID** and click **OK** to open the Analysis Method Editor with the General tab selected.
Enter the settings shown in the figures on the following pages.
4. After you enter settings in all tabs, click **Save**.

General tab settings



In the Name field, either type the name as shown for consistency with files supplied with other AmpF Φ STR™ kits, or enter a name of your choosing. The Description and Instrument fields are optional.

Allele tab settings



- In the Bin Set field, select the **NGM_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.
- The “Use marker-specific stutter ratio if available” check box is selected by default. Consequently, the software applies the stutter ratio filters supplied in the NGM_panel_v2 file. GeneMapper™ ID Software v3.2.1 specifies locus-specific filter ratios for minus stutters, but not for plus stutters, in the panel file. However, validation studies with the NGM™ Kit show that the trinucleotide repeat D22S1045 locus produces a relatively large amount of plus stutter compared to tetranucleotide repeat loci. The relatively large amount of stutter may cause the stutter peak to be labeled during routine analysis.
- The plus stutter at the D22S1045 locus can be filtered by assigning a global plus stutter filter for trinucleotide repeat loci in the Analysis Parameter file. Because D22S1045 is the only trinucleotide repeat locus in the NGM™ Kit, this stutter filter setting is applied only to plus stutter peaks at the D22S1045 locus. The settings shown above resulted in little or no labeling of D22S1045 plus stutter peaks during our validation studies. Perform internal validation studies to determine the settings to use in your laboratory.

Peak Detector tab settings

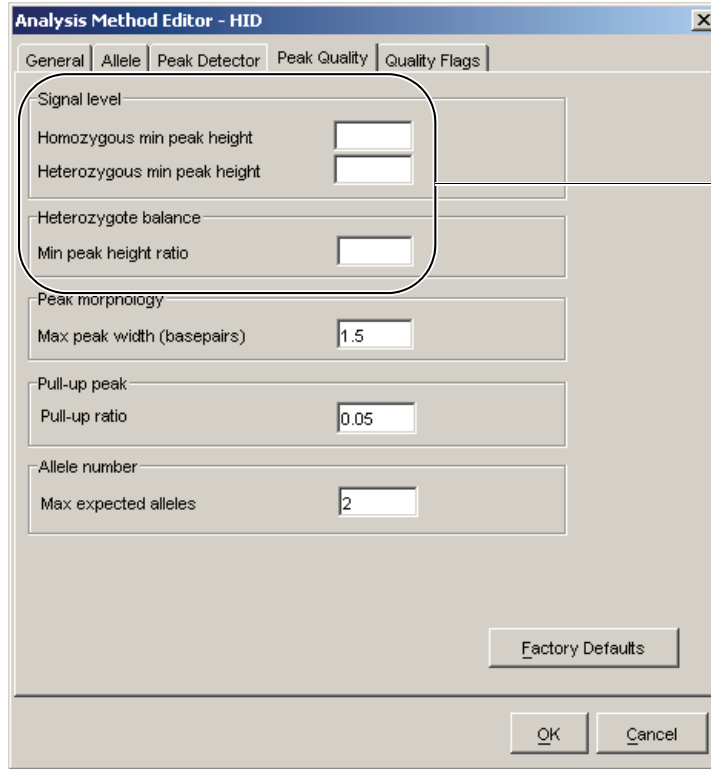
The screenshot shows the 'Analysis Method Editor - HID' window with the 'Peak Detector' tab selected. The 'Peak Amplitude Thresholds' section is circled in red, and a callout box points to it with the text 'Perform internal validation studies to determine settings'. The 'Peak Amplitude Thresholds' section includes fields for B, R, G, O, and Y. Other settings include 'Peak Detection Algorithm' set to 'Advanced', 'Analysis' set to 'Full Range', 'Sizing' set to 'All Sizes', 'Smoothing' set to 'Light', 'Baseline Window' set to '51 pts', and 'Size Calling Method' set to '3rd Order Least Squares'.

IMPORTANT! Perform the appropriate internal validation studies to determine the peak amplitude thresholds for interpretation of NGM™ 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 NGM™ Kit has been validated using the 3rd Order Least Squares sizing method in combination with the GeneScan™ 500 LIZ™ Size Standard. If using the GeneScan™ 600 LIZ™ Size Standard v2.0, select the Local Southern Method. Select alternative sizing methods only after you perform the appropriate internal validation studies.

Peak Quality tab settings



Perform internal validation studies to determine settings

IMPORTANT! Perform the appropriate internal validation studies to determine the heterozygous and homozygous minimum peak height thresholds and the minimum peak height ratio threshold that allow for reliable interpretation of NGM™ Kit data.

Quality Flags tab settings

Analysis Method Editor - HID

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

Quality weights are between 0 and 1.

Quality Flag Settings

Spectral Pull-up	0.8	Control Concordance	1.0
Broad Peak	0.8	Low Peak Height	0.3
Out of Bin Allele	0.8	Off-scale	0.8
Overlap	0.8	Peak Height Ratio	0.3

P/QV Thresholds

	Pass Range:	Low Quality Range:
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

Factory Defaults

OK Cancel

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 to use in your laboratory.

Create a size standard

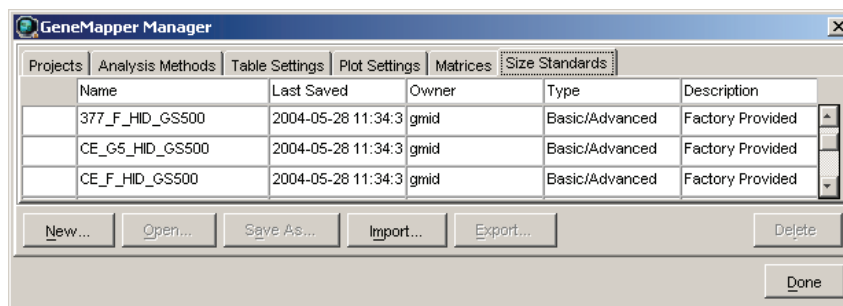
The size standards for the NGM™ Kit uses the following size standard peaks in their definitions:

GeneScan™ 500 LIZ™ Size Standard	GeneScan™ 600 LIZ™ Size Standard v2.0
75, 100, 139, 150, 160, 200, 300, 350, 400, and 450	80, 100, 114, 120, 140, 160, 180, 200, 214, 220, 240, 250, 260, 280, 300, 314, 320, 340, 360, 380, 400, 414, 420, 440 and 460

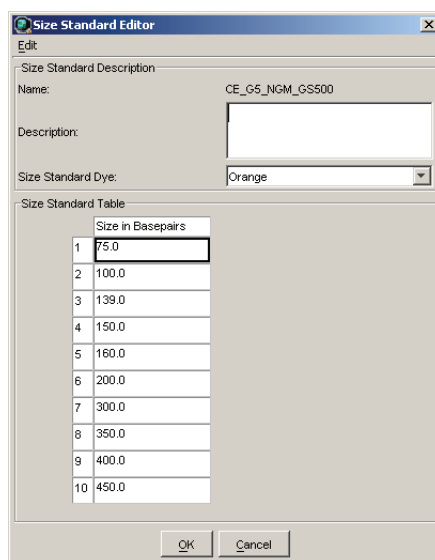
Note: The 250-nt and the 340-nt peak are not included in the size standard definition. These peaks can be used as an indicator of precision within a run.

Use the following procedure to create the size standard for the NGM™ Kit.

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



2. Select the **Size Standards** tab, then click **New**.
3. Enter a name as shown below or enter a name of your choosing. In the Size Standard Dye field, select **Orange**. In the Size Standard Table, enter the sizes specified in on page 42. The example below is for the GeneScan™ 500 LIZ™ Size Standard.




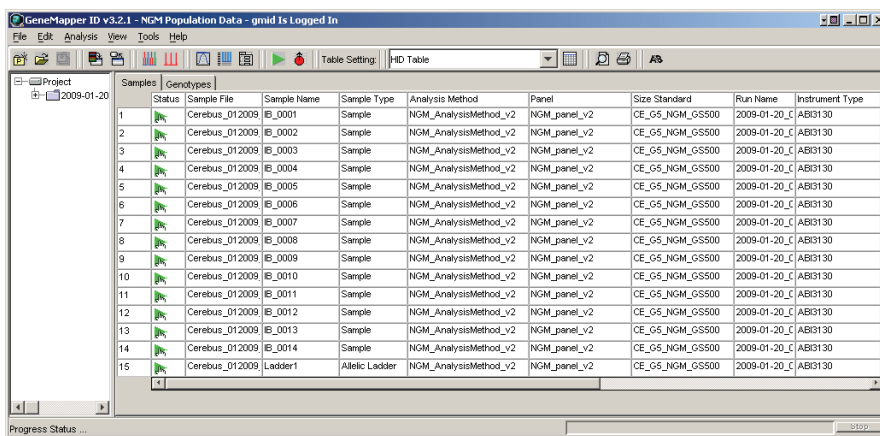
Analyze and edit sample files with GeneMapper™ ID Software

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. The names of the settings shown are the names suggested in the sections above.

Parameter	Settings
Sample Type	Select the sample type.
Analysis Method	NGM_AnalysisMethod_v2 (or the name of the analysis method you created)
Panel	NGM_panel_v2
Size Standard	CE_G5_NGM_GS500 (or the name of the size standard you created)

For more information about how the Size Caller works, refer to the ABI Prism™ GeneScan™ Analysis Software for the Windows NT™ Operating System Overview of the Analysis Parameters and Size Caller User Bulletin (Part no. 4335617). For additional information about size standards, refer to the GeneMapper™ ID Software Version 3.1 Human Identification Analysis User Guide (Part no. 4338775).

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 both:
 - A completion bar extending to the right with the percentage completed 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.



The screenshot shows the GeneMapper ID v3.2.1 software interface. The main window displays a table with columns: Status, Sample File, Sample Name, Sample Type, Analysis Method, Panel, Size Standard, Run Name, and Instrument Type. The table contains 15 rows of data, with the first row highlighted in green. The status bar at the bottom shows 'Progress Status ...' and a 'stop' button.

Status	Sample File	Sample Name	Sample Type	Analysis Method	Panel	Size Standard	Run Name	Instrument Type
1	Cerebus_012009_E_0001		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
2	Cerebus_012009_E_0002		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
3	Cerebus_012009_E_0003		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
4	Cerebus_012009_E_0004		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
5	Cerebus_012009_E_0005		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
6	Cerebus_012009_E_0006		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
7	Cerebus_012009_E_0007		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
8	Cerebus_012009_E_0008		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
9	Cerebus_012009_E_0009		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
10	Cerebus_012009_E_0010		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
11	Cerebus_012009_E_0011		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
12	Cerebus_012009_E_0012		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
13	Cerebus_012009_E_0013		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
14	Cerebus_012009_E_0014		Sample	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130
15	Cerebus_012009_Ladder1		Allelic Ladder	NGM_AnalysisMethod_v2	NGM_panel_v2	CE_G5_NGM_GS500	2009-01-20_C	ABI3130

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

Note: .hid files can only be analyzed using GeneMapper™ ID-X Software v1.2 or higher

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. Perform the appropriate internal validation studies if you want to use multiple ladder samples in an analysis.
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 ± 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, verify the sample result according to your laboratory protocol.

Set up GeneMapper™ ID-X 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.appliedbiosystems.com.

Overview To analyze sample (.fsa) files using GeneMapper™ ID-X Software v1.0.1 or higher for the first time:

1. Import panels, bins, and marker stutter into the Panel Manager, as explained in “Import panels, bins, and marker stutter” below.
2. Create an analysis method, as explained in “Create an analysis method” on page 50.
3. Create a size standard, as explained in “Create a size standard” on page 55.
4. Define custom views of analysis tables.
Refer to the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Part no. 4375574) for more information.
5. Define custom views of plots.
Refer to the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Part no. 4375574) for more information.

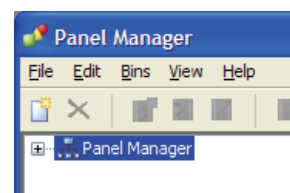
Import panels, bins, and marker stutter

To import the NGM™ Kit panels, bin sets, and marker stutter from the Applied Biosystems web site into the GeneMapper™ ID-X Software database:

1. Download and open the file containing panels, bins, and marker stutter:
 - a. From the Support menu of www.appliedbiosystems.com, select **Support ▶ Software Downloads, Patches & Updates ▶ GeneMapper™ ID-X Software ▶ Updates & Patches**, and download the file **NGM Analysis Files GMIDX**.
 - b. Unzip the file.
2. Start the GeneMapper™ ID-X Software, then log in with the appropriate user name and password.

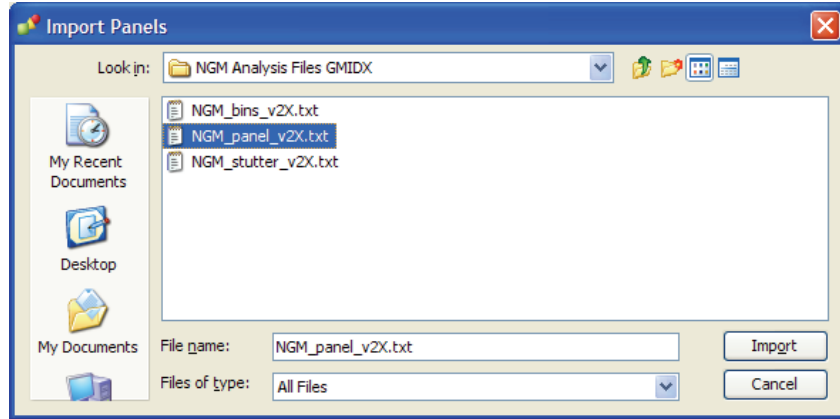
IMPORTANT! For logon instructions, refer to the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Part no. 4375574).

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 **NGM Analysis Files GMIDX** folder that you unzipped in step 1 of this section.

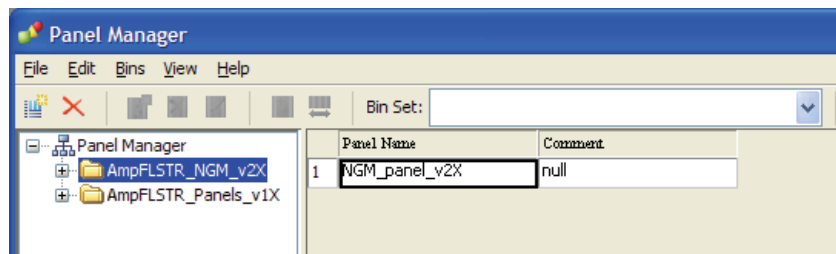


5. Select **NGM_panel_v2X**, then click **Import**.

Note: Importing this file creates a new folder in the navigation pane of the Panel Manager “AmpFLSTR_NGM_v2X”. This folder contains the panel and associated markers.

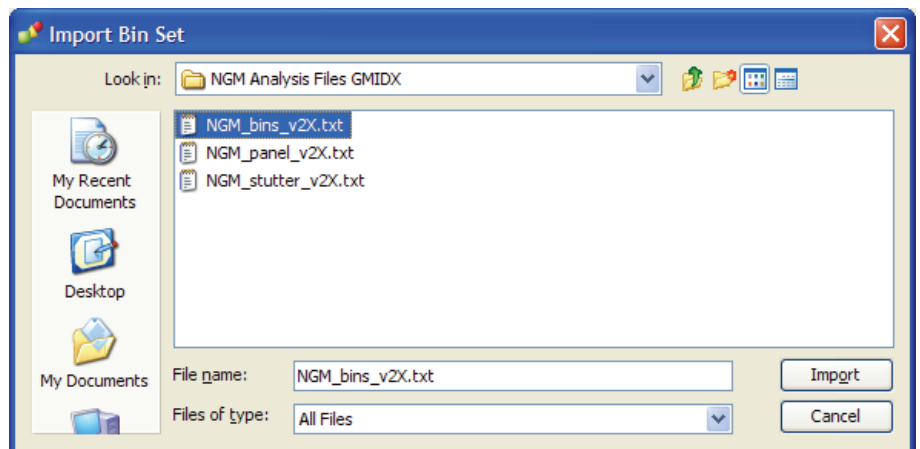


6. Import **NGM_bins_v2X.txt**:
 - a. Select the **AmpFLSTR_NGM_v2X** folder in the navigation pane.



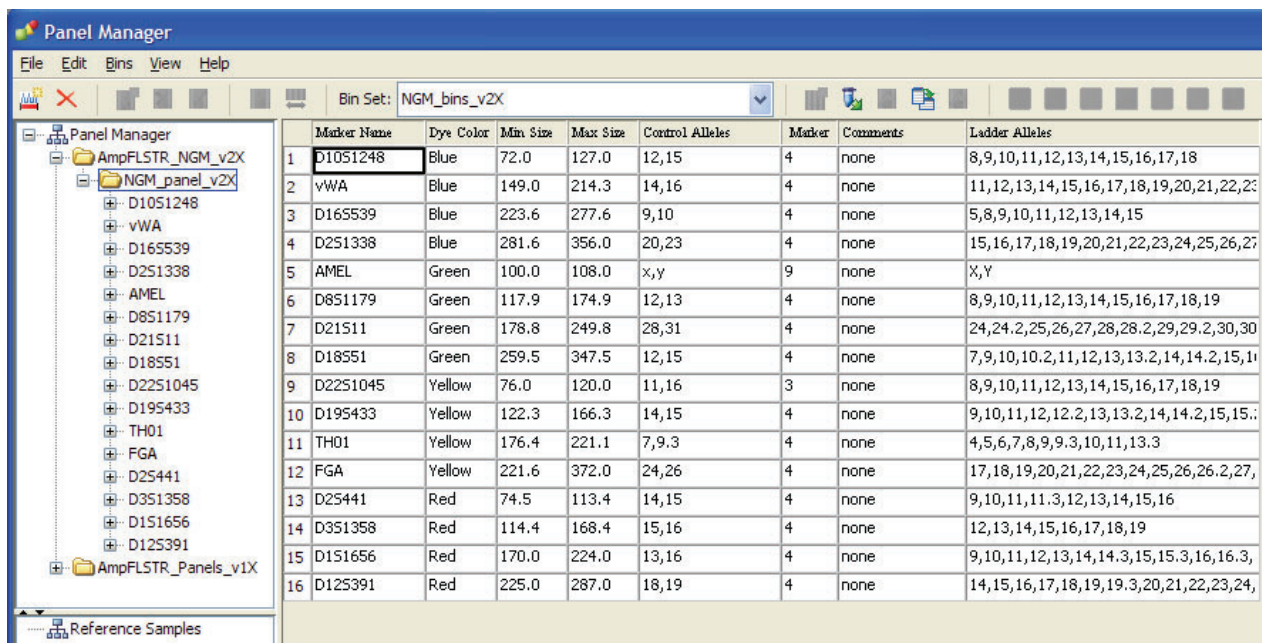
- b. Select **File** ► **Import Bin Set** to open the Import Bin Set dialog box.
 - c. Navigate to, then open the **NGM Analysis Files GMIDX** folder.
 - d. Select **NGM_bins_v2X.txt**, then click **Import**.

Note: Importing this file associates the bin set with the panels in the **NGM_panel_v2X** folder.



7. View the imported panels in the navigation pane:

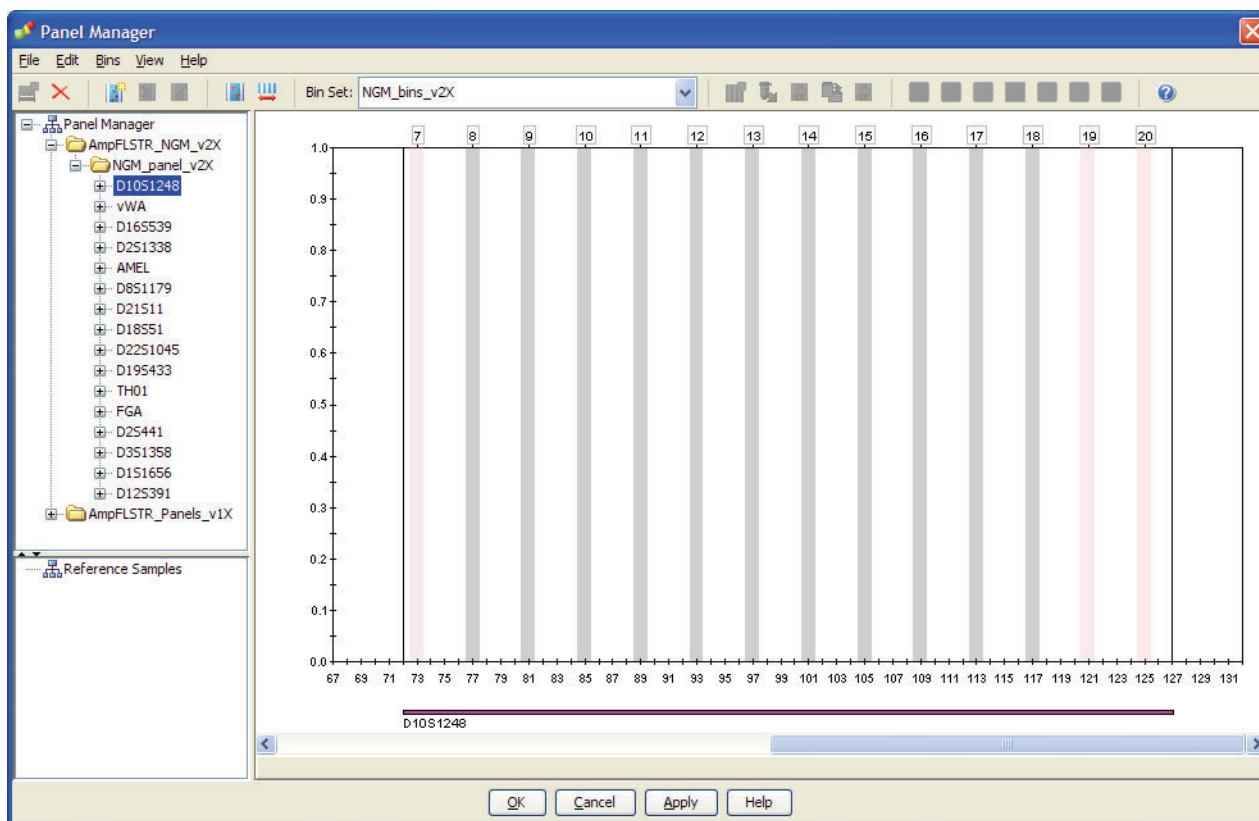
- a. Double-click the **AmpFLSTR_NGM_v2X** folder.
- b. Double-click the **NGM_panel_v2X** folder to display the panel information in the right pane and the markers below it.



The screenshot shows the Panel Manager software interface. The left pane displays a tree view with folders: AmpFLSTR_NGM_v2X, NGM_panel_v2X, and AmpFLSTR_Panels_v1X. The right pane shows a table of markers with the following columns: Marker Name, Dye Color, Min. Size, Max. Size, Control Alleles, Marker, Comments, and Ladder Alleles. The marker D10S1248 is highlighted in the table.

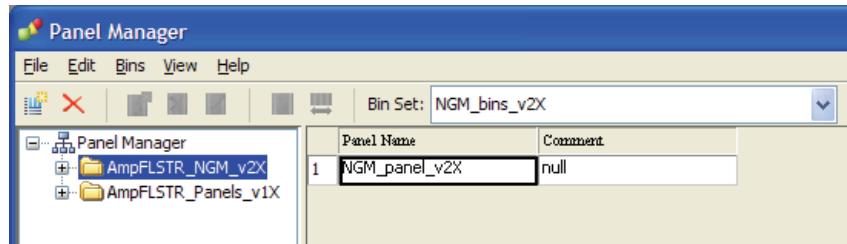
Marker Name	Dye Color	Min. Size	Max. Size	Control Alleles	Marker	Comments	Ladder Alleles
1 D10S1248	Blue	72.0	127.0	12,15	4	none	8,9,10,11,12,13,14,15,16,17,18
2 vWA	Blue	149.0	214.3	14,16	4	none	11,12,13,14,15,16,17,18,19,20,21,22,23
3 D16S539	Blue	223.6	277.6	9,10	4	none	5,8,9,10,11,12,13,14,15
4 D2S1338	Blue	281.6	356.0	20,23	4	none	15,16,17,18,19,20,21,22,23,24,25,26,27
5 AMEL	Green	100.0	108.0	x,y	9	none	X,Y
6 D8S1179	Green	117.9	174.9	12,13	4	none	8,9,10,11,12,13,14,15,16,17,18,19
7 D21S11	Green	178.8	249.8	28,31	4	none	24,24.2,25,26,27,28,28.2,29,29.2,30,30.2
8 D18S51	Green	259.5	347.5	12,15	4	none	7,9,10,10.2,11,12,13,13.2,14,14.2,15,15.2
9 D22S1045	Yellow	76.0	120.0	11,16	3	none	8,9,10,11,12,13,14,15,16,17,18,19
10 D19S433	Yellow	122.3	166.3	14,15	4	none	9,10,11,12,12.2,13,13.2,14,14.2,15,15.2
11 TH01	Yellow	176.4	221.1	7,9,3	4	none	4,5,6,7,8,9,9.3,10,11,13.3
12 FGA	Yellow	221.6	372.0	24,26	4	none	17,18,19,20,21,22,23,24,25,26,26.2,27,27.2
13 D2S441	Red	74.5	113.4	14,15	4	none	9,10,11,11.3,12,13,14,15,16
14 D3S1358	Red	114.4	168.4	15,16	4	none	12,13,14,15,16,17,18,19
15 D1S1656	Red	170.0	224.0	13,16	4	none	9,10,11,12,13,14,14.3,15,15.3,16,16.3,17
16 D12S391	Red	225.0	287.0	18,19	4	none	14,15,16,17,18,19,19.3,20,21,22,23,24,25

8. Select **D10S1248** to display the Bin view for the marker in the right pane.



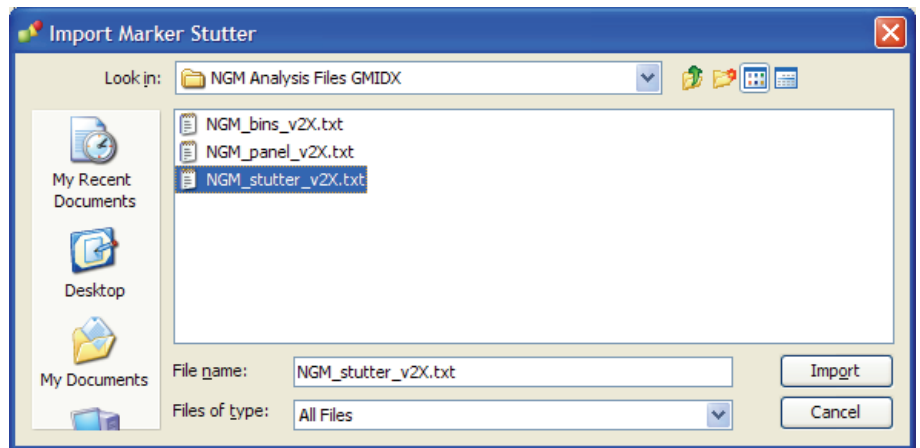
9. Import **NGM_stutter_v2X.txt**:

- a. Select the **AmpFLSTR_NGM_v2** folder in the navigation panel.



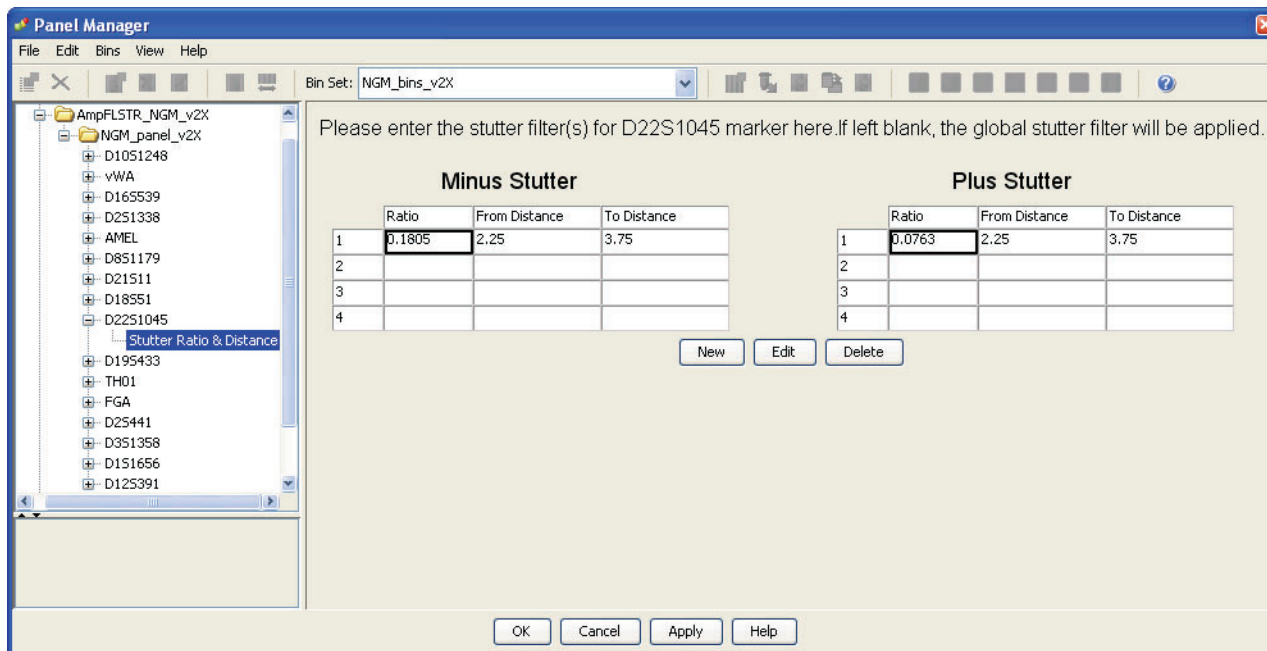
- b. Select File ► **Import Marker Stutter** to open the Import Marker Stutter dialog box.
- c. Navigate to, then open the **NGM Analysis Files GMIDX** folder.
- d. Select **NGM_stutter_v2X.txt**, then click **Import**.

Note: Importing this file associates the marker stutter ratio with the bin set in the NGM_bins_v2X folder.



10. View the imported marker stutters in the navigation pane:
 - a. Select the **NGM_panel_v2X** folder to display its list of markers in the right pane.
 - b. Double-click the **NGM_panel_v2X** folder to display its list of markers below it.
 - c. Double-click **D22S1045** to display the Stutter Ratio & Distance view for the marker in the right pane.

Because D22S1045 has a trinucleotide repeat unit, it produces a higher level of plus stutter than tetranucleotide markers, and so requires the use of a plus stutter filter. The settings for the D22S1045 plus stutter filter can be seen in the table in the right pane. Other markers may not require a plus stutter filter, in which case the settings for plus stutter are left blank.



11. Click **Apply**, then **OK** to add the NGM™ Kit panels, bin sets, and marker stutter to the GeneMapper™ ID-X Software database.

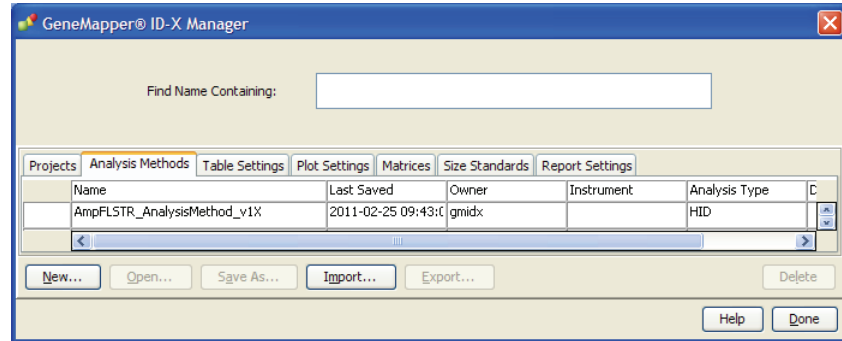
IMPORTANT! If you close the Panel Manager without clicking **Apply**, the panels, bin sets, and marker stutter will not be imported into the GeneMapper™ ID-X Software database.

Create an analysis method

Use the following procedure to create an analysis method for the NGM™ 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 Software 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.



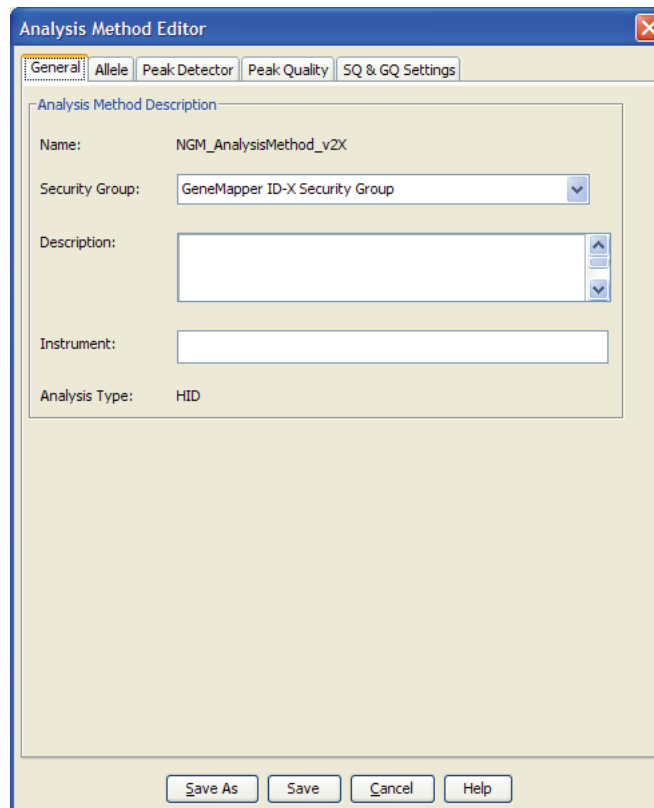
2. Select the **Analysis Methods** tab, then click **New** to open the Analysis Method Editor with the **General** tab selected.

3. Enter the settings shown in the figures on the following pages.

Note: The Analysis Method Editor closes when you save your settings (see step 4 on page 51). To complete this step quickly, do not save the analysis method until you finish entering settings in all of the tabs.

4. After you enter the settings on all tabs, click **Save**.

General tab settings



In the Name field, either type the name as shown for consistency with files supplied with other AmpFISTR™ kits or enter a name of your choosing. In the Security Group field, select the Security Group appropriate to your software configuration from the drop-down list. The Description and Instrument fields are optional.

Allele tab settings

Analysis Method Editor

General Allele Peak Detector Peak Quality SQ & GQ Settings

Bin Set: NGM_bins_v2X

Use marker-specific stutter ratio and distance if available

Marker Repeat Type:		Tri	Tetra	Penta	Hexa
Global 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
Global Minus Stutter Ratio		0.0	0.0	0.0	0.0
Global Minus Stutter Distance	From	2.25	3.25	0.0	0.0
	To	3.75	4.75	0.0	0.0
Global Plus Stutter Ratio		0.0	0.0	0.0	0.0
Global Plus Stutter Distance	From	0.0	0.0	0.0	0.0
	To	0.0	0.0	0.0	0.0

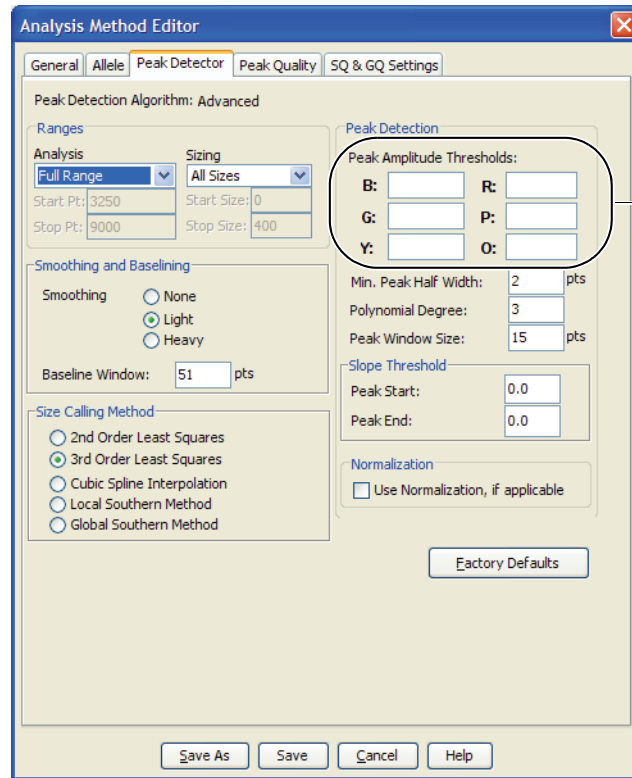
Amelogenin Cutoff: 0.0

Range Filter... Factory Defaults

Save As Save Cancel Help

- In the Bin Set field, select the **NGM_bins_v2X** bin set imported previously and configure the stutter distance parameters as shown.
- GeneMapper™ ID-X Software 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.
- The “Use marker-specific stutter ratio if applicable” check box is selected by default. When this box is checked, the software applies the stutter ratio filters in the NGM_stutter_v2X.txt file.

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 NGM™ 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 NGM™ Kit has been validated using the 3rd Order Least Squares sizing method with the GeneScan™ 500 LIZ™ Size Standard. If you use GeneScan™ 600 LIZ™ Size Standard v2.0, select the Local Southern Method. Select alternative sizing methods only after you perform the appropriate internal validation studies.
- **Normalization (v1.2 or higher)** – For use with 3500 data. Perform internal validation studies to determine whether to use the Normalization feature for analysis of NGM™ 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

Spike Detection

Cut-off Value

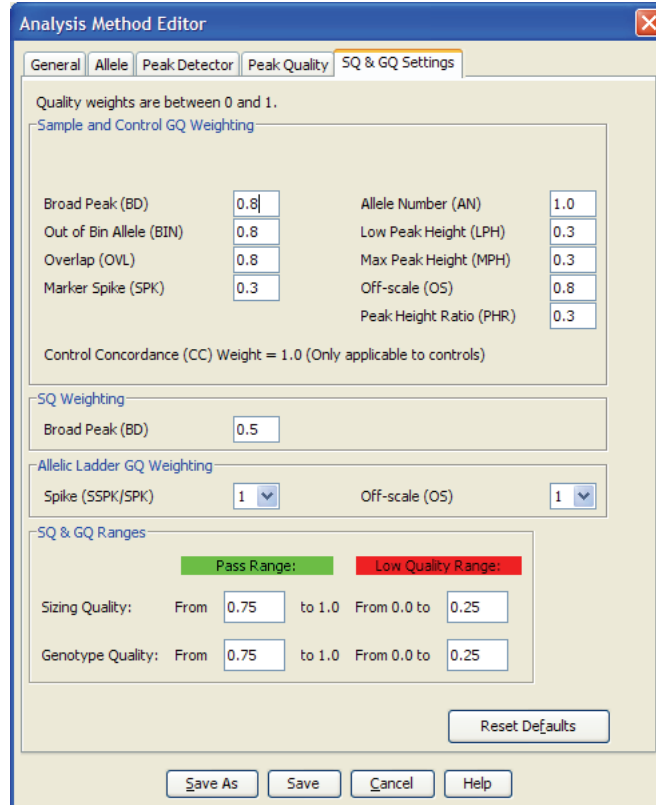
Factory Defaults

Save As Save Cancel Help

Perform internal validation studies to determine settings

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

SQ & GQ tab settings



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 a size standard

The size standards for the NGM™ Kit uses the following size standard peaks in their definitions:

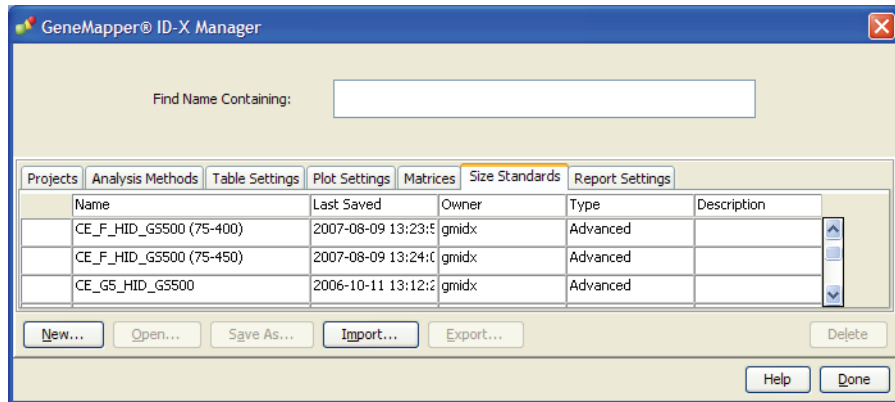
GeneScan™ 500 LIZ™ Size Standard	GeneScan™ 600 LIZ™ Size Standard v2.0
75, 100, 139, 150, 160, 200, 300, 350, 400, and 450	80, 100, 114, 120, 140, 160, 180, 200, 214, 220, 240, 250, 260, 280, 300, 314, 320, 340, 360, 380, 400, 414, 420, 440 and 460

Note: The 250-nt and the 340-nt peaks are not included in the size standard definition. These peaks can be used as an indicator of precision within a run.

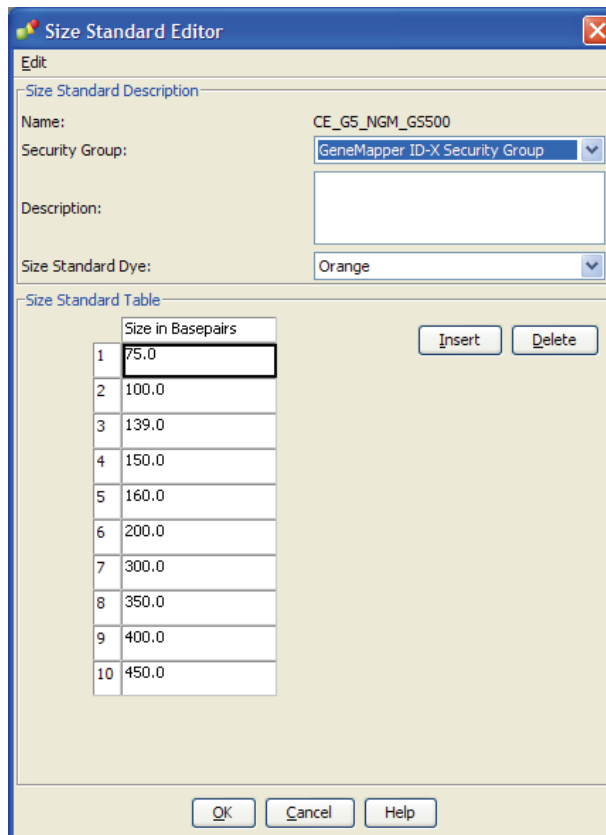
Use the following procedure to create the size standard for the NGM™ Kit.

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

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



3. Complete the Name field as shown below or with a name of your choosing. In the Security Group field, select the Security Group appropriate to your software configuration from the drop-down list. In the Size Standard Dye field, select **Orange**. In the Size Standard Table, enter the sizes specified on page 55. The example below is for the GeneScan™ 500 LIZ™ Size Standard.




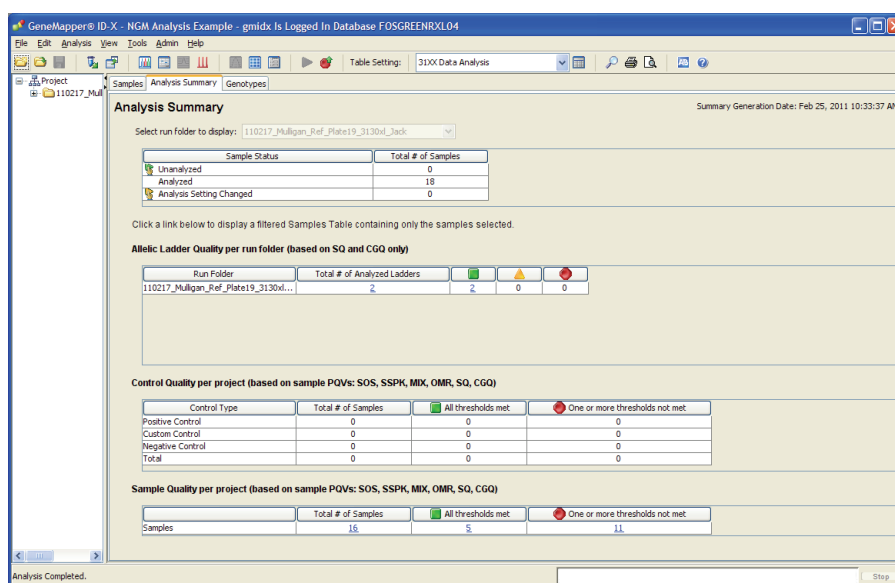
Analyze and edit sample files with GeneMapper™ ID-X Software

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. The names of the settings shown are the names suggested in the sections above. If you named the settings differently, select the names you specified.

Parameter	Settings
Sample Type	Select the sample type.
Analysis Method	NGM_AnalysisMethod_v2X (or the name of the analysis method you created)
Panel	NGM_panel_v2X
Size Standard	CE_G5_NGM_GS500 (or the name of the size standard you created)

For more information about how the Size Caller works, or about size standards, refer to the *GeneMapper™ ID-X Software v1.2 Reference Guide* (Part no. 4426481A).

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 completed 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 upon completion of the analysis.
- The figure below shows the 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 about any of these tasks, refer to:

- For quick set-up instructions, refer to the *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Part no. 4375574).
- For details about GeneMapper™ ID-X Software features, allele filters, peak detection algorithms, and project editing, refer to:
 - *GeneMapper™ ID-X Software Version 1.0 Getting Started Guide* (Part no. 4375574)
 - *GeneMapper™ ID-X Software Version 1.0 Quick Reference Guide* (Part no. 4375670)
 - *GeneMapper™ ID-X Software Version 1.0 Reference Guide* (Part no. 4375671)
 - *GeneMapper™ ID-X Software Version 1.1 (Mixture Analysis Tool) Getting Started Guide* (Part no. 4396773)
 - *GeneMapper™ ID-X Software Version 1.1 (Mixture Analysis Tool) Quick Reference Guide* (Part no. 4402094)
 - *GeneMapper™ ID-X Software Version 1.2 Quick Reference Guide* (Part no. 4426482)
 - *GeneMapper™ ID-X Software Version 1.2 Reference Guide* (Part no. 4426481)

■ Overview	59
■ Developmental validation.....	60
■ Accuracy, precision, and reproducibility	63
■ Extra peaks in the electropherogram.....	72
■ Characterization of loci	79
■ Species specificity.....	81
■ Sensitivity	82
■ Stability	84
■ Mixture studies	88
■ Population data.....	91
■ New Primers added to the NGM™ Kit	91
■ Mutation rate.....	93
■ Probability of identity	94
■ Probability of paternity exclusion	104

Overview

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 that 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

We performed experiments to evaluate the performance of the NGM™ Kit. The experiments were performed according to the revised guidelines from the Scientific Working Group on DNA Analysis Methods (SWGDM, July 10, 2003). Based on these guidelines, we conducted experiments that comply with guidelines 1.0 and 2.0 and its associated subsections. This DNA methodology is not novel. (Moretti *et al.*, 2001; Frank *et al.*, 2001; Wallin *et al.*, 2002; and Holt *et al.*, 2000).

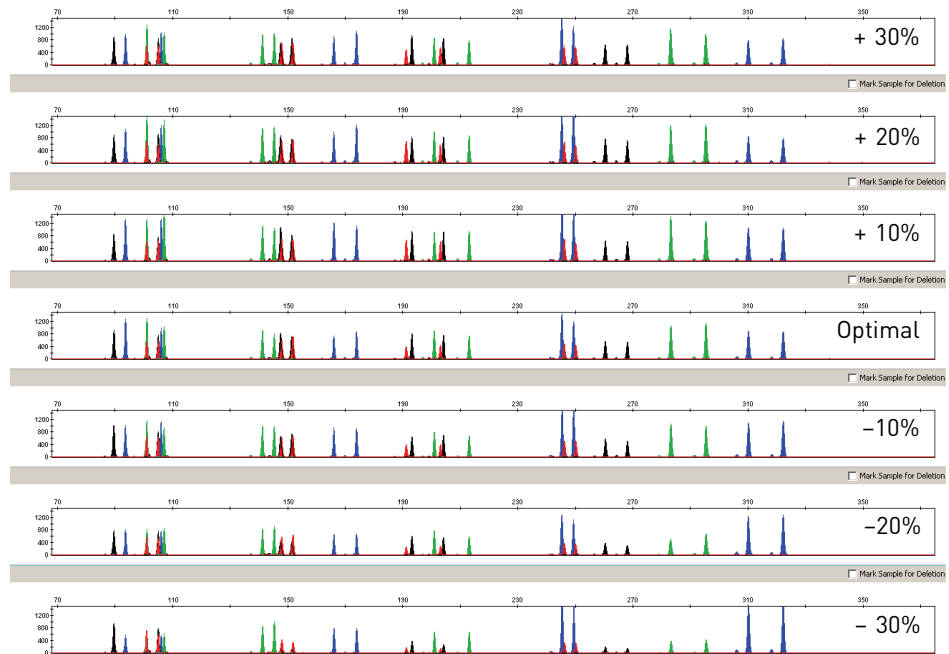
This chapter discusses many of the experiments performed by us and provides examples of results obtained. We chose conditions that produced optimum PCR product yield and that met reproducible performance standards. It is our opinion that while these experiments are not exhaustive, they are appropriate for a manufacturer of STR kits intended for forensic and/or parentage testing use.

IMPORTANT! Perform internal validation studies before using the NGM™ Kit.

Developmental validation

- SWGAM guideline 1.2.1** “Developmental validation is the demonstration of the accuracy, precision, and reproducibility of a procedure by the manufacturer, technical organization, academic institution, government laboratory, or other party.” (SWGAM, July 2003)
- SWGAM guideline 2.10.1** “The reaction conditions needed to provide the required degree of specificity and robustness must be determined. These include thermocycling parameters, the concentration of primers, magnesium chloride, DNA polymerase, and other critical reagents.” (SWGAM, July 2003)
- PCR components** We examined the concentration of each component of the NGM™ Kit and established that the concentration of each component was within the range where data indicated that the amplification met the required performance criteria for specificity, sensitivity, and reproducibility. For example, 1.0 ng of Control DNA 007 was amplified in the presence of varying concentrations of magnesium chloride, and the results were analyzed on an Applied Biosystems™ 3130xl Genetic Analyzer. Results are shown in Figure 5. The performance of the multiplex is most robust within ±20% of the optimal magnesium chloride concentration.

Figure 5 1.0 ng of control DNA 007 amplified for 29 cycles with the NGM™ Kit in the presence of varying concentrations of magnesium chloride and analyzed on an Applied Biosystems™ 3130xl Genetic Analyzer



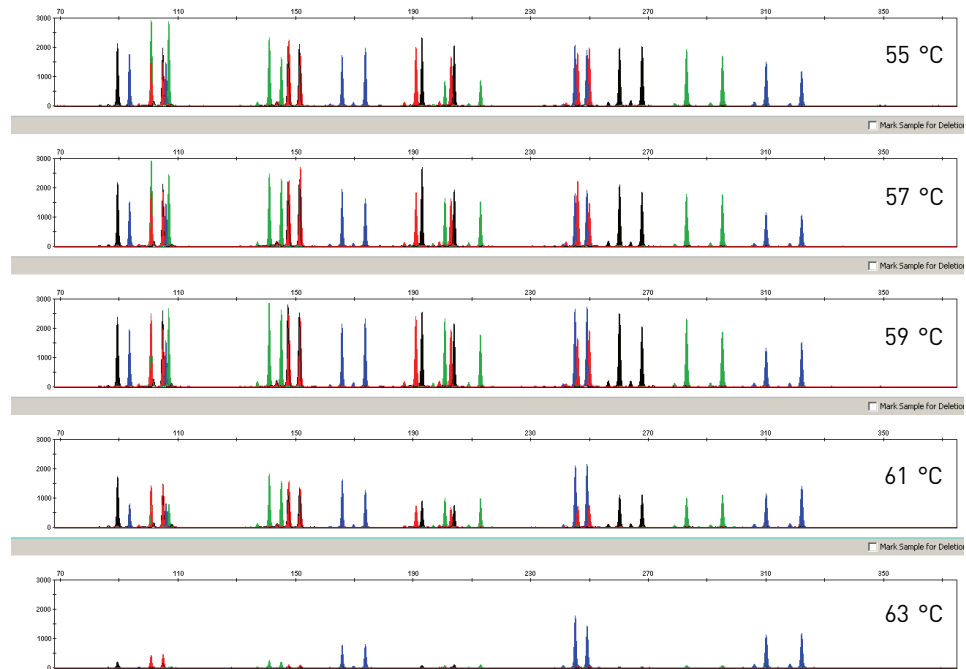
Thermal cycler parameters

Thermal cycling parameters were optimized using a Design of Experiments (DOE) approach that attempts to identify the combination of temperatures and hold times that produce the best assay performance. Optimal assay performance was determined through evaluation of assay sensitivity, peak-height balance and resistance to PCR inhibitors.

For example, annealing temperatures of 55, 57, 59, 61, and 63°C were tested using a Silver 96-Well GeneAmp™ PCR System 9700 (Figure 6). The PCR products were analyzed using an Applied Biosystems™ 3130xl Genetic Analyzer.

Of the tested annealing temperatures, 55 to 61°C produced robust profiles. At 63°C the yield of the majority of loci was significantly reduced. The optimal combination of specificity, sensitivity, and resistance to PCR inhibition was observed at 59°C. Thermal cycler temperature is critical to assay performance; therefore, routine, regularly scheduled thermal cycler calibration is strongly recommended.

Figure 6 Electropherograms obtained from amplification of 1.0 ng of control DNA 007 for 29 cycles at annealing temperatures of 55, 57, 59, 61, and 63°C, analyzed on an Applied Biosystems™ 3130xl Genetic Analyzer, (Y-axis scale 0 to 3,000 RFU)

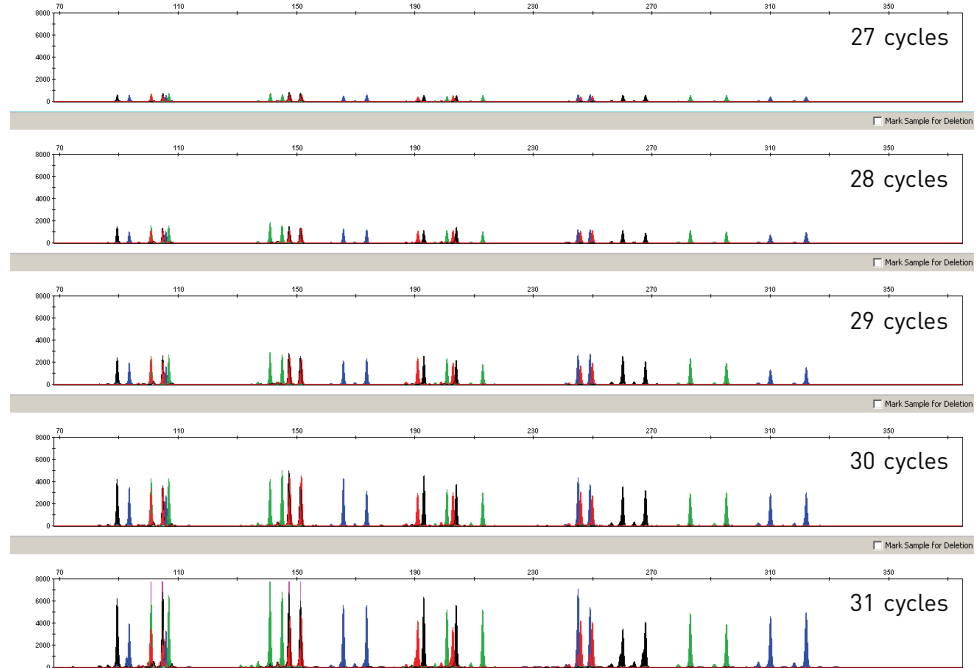


PCR cycle number

NGM™ Kit reactions were amplified for 27, 28, 29, 30, and 31 cycles on the Silver 96-Well GeneAmp™ PCR System 9700 using 1.0 ng of each of three DNA samples. As expected, the amount of PCR product increased with the number of cycles. A full profile was generated for all numbers of thermal cycles (27-31) and off-scale data were collected for several allele peaks at 30 and 31 cycles (Figure 7).

The NGM™ Kit was optimized for use with both 29 or 30 cycles and validated on the Applied Biosystems™ 3130xl Genetic Analyzer. The optimal DNA input amount per PCR was found to be 1.0 ng and 0.5 ng, respectively, for 29- and 30-cycle amplifications. None of the cycle numbers tested produced nonspecific peaks.

Figure 7 Representative NGM™ Kit profiles obtained from amplification of 1.0 ng DNA template using 27, 28, 29, 30, and 31 cycles, analyzed on an Applied Biosystems™ 3130xl Genetic Analyzer, (Y-axis scale 0 to 8,000 RFU)



Accuracy, precision, and reproducibility

SWGDM guideline 2.9

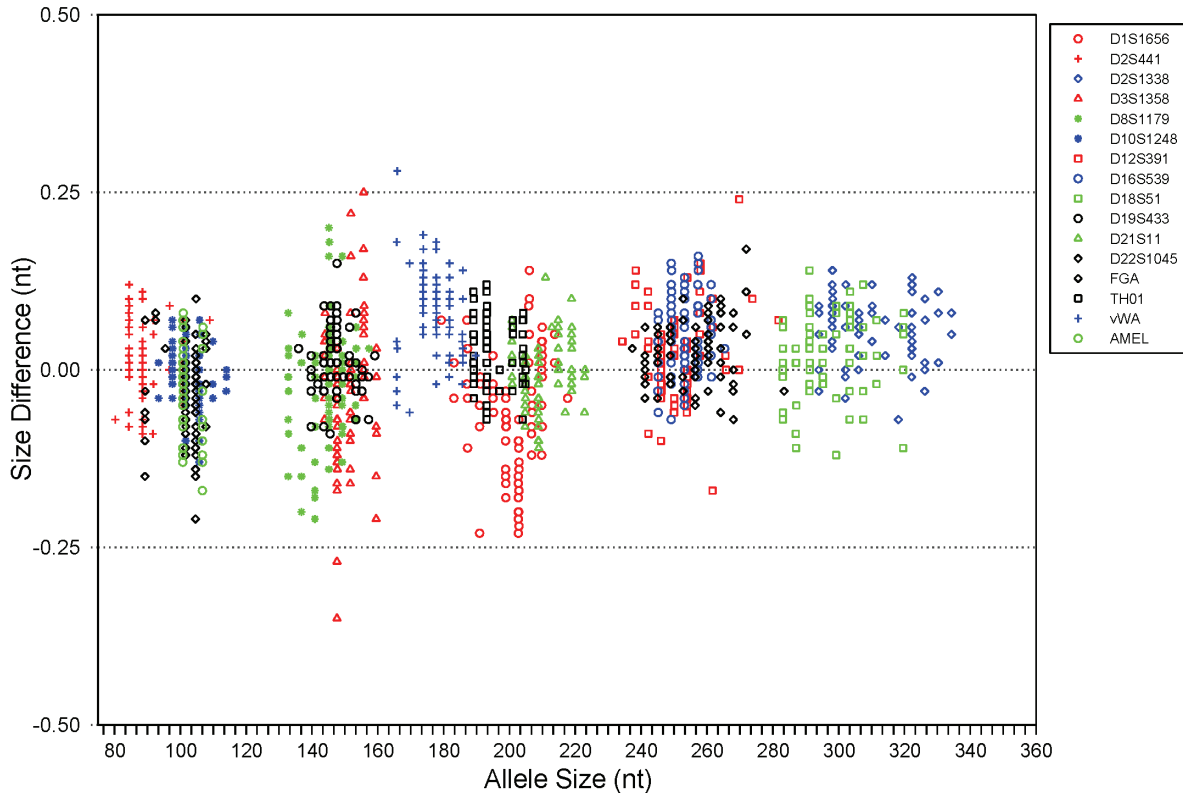
“The extent to which a given set of measurements of the same sample agree with their mean and the extent to which these measurements match the actual values being measured should be determined.” (SWGDM, July 2003)

Accuracy

Laser-induced fluorescence detection of length polymorphism at short tandem repeat loci is not a novel methodology (Holt *et al.*, 2000; and Wallin *et al.*, 2002). However, accuracy and reproducibility of NGM™ Kit profiles have been determined from various sample types. Figure 8 shows the size differences that are typically observed between sample alleles and allelic ladder alleles on the Applied Biosystems™ 3130xl Genetic Analyzer with POP-4™ polymer. The x-axis in Figure 8 represents the nominal nucleotide sizes for the AmpFSTR™ NGM™ Allelic Ladder. The dashed lines parallel to the x-axis represent the ± 0.25 -nt windows. The y-axis represents the deviation of each sample allele size from the corresponding allelic ladder allele size. All sample alleles are within ± 0.5 nt from a corresponding allele in the allelic ladder.

Figure 8 Allele Size vs. Allelic Ladder Sizing for 42 samples analyzed on an Applied Biosystems™ 3130xl Genetic Analyzer.

Size and ladder sizing for the NGM™ Kit were calculated using the GeneScan™ 500 LIZ™ Size Standard using the 3rd Order Least Squares Method



Precision and size windows

Sizing precision enables the determination of accurate and reliable genotypes. Sizing precision was measured on an Applied Biosystems™ 3130xl Genetic Analyzer. The recommended method for genotyping is to employ a ± 0.5 -nt “window” around the size obtained for each allele in the AmpF ℓ STR™ NGM™ Allelic Ladder. A ± 0.5 -nt window allows for the detection and correct assignment of alleles. Any sample allele that sizes outside the specified window could be:

- An “off-ladder” allele, that is, an allele of a size that is not represented in the AmpF ℓ STR™ NGM™ Allelic Ladder

or

- 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 on a capillary instrument or in several lanes of one gel.

Table 7 on page 86 shows typical precision results obtained from five runs (16 capillaries/run) of the AmpF ℓ STR™ NGM™ Allelic Ladder on an Applied Biosystems™ 3130xl Genetic Analyzer (36-cm capillary and POP-4™ polymer), using the GeneScan™ 500 LIZ™ Size Standard. The results were obtained within a set of injections on a single capillary array.

Sample alleles may occasionally size outside of the ± 0.5 -nt window for a respective allelic ladder allele because of measurement error. The frequency of such an occurrence is lowest in detection systems having the smallest standard deviations in sizing. Figure 8 on page 64 illustrates the tight clustering of allele sizes obtained on the Applied Biosystems™ 3130xl Genetic Analyzer, where the standard deviation in sizing is typically less than 0.15 nt. The instance of a sample allele sizing outside the ± 0.5 -nt window because of measurement error is relatively rare when the standard deviation in sizing is approximately 0.15 nt or less (Smith, 1995).

For sample alleles that do not size within a ± 0.5 -nt 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. Repeat analysis, when necessary, provides an added level of confidence in the final allele assignment.

GeneMapper™ ID Software and GeneMapper™ ID-X Software automatically flag sample alleles that do not size within the prescribed window around an allelic ladder allele by labelling the allele as OL (off-ladder).

Maximum sizing precision is obtained within the same set of capillary injections. Cross-platform sizing differences occur due to a number of factors including type and concentration of polymer, run temperature, and electrophoresis conditions. Variations in sizing can also occur between runs on the same instrument and between runs on different instruments of the same platform type because of these factors.

We strongly recommend that the allele sizes be compared to the sizes obtained for known alleles in the AmpF Φ STR™ NGM™ Allelic Ladder from the same run and then be converted to genotypes (as described in “Before you start” on page 34 and on page 45). See Table 4 for the results of five runs of the AmpF Φ STR™ NGM™ Allelic Ladder on an Applied Biosystems™ 3130xl Genetic Analyzer. For more information on precision and genotyping, see Lazaruk *et al.*, 1998 and Mansfield *et al.*, 1998.

In Table 4, the mean sizes for all the alleles in each run (16 capillaries) were calculated. The mean range shown in the table represents the lowest- and highest-mean size values obtained across all five runs. Similarly, the standard deviation for the allele sizing was calculated for all the alleles in each run. The standard deviation range shown in Table 4 represents the lowest and highest standard deviation values obtained across all five runs.

Table 4 Precision results of five runs (16 capillaries/run) of the AmpF Φ STR™ NGM™ Allelic Ladder

Allele	Mean	Standard Dev.
Amelogenin		
X	100.54–100.68	0.052–0.064
Y	106.21–106.34	0.045–0.061
D10S1248		
8	76.64–76.67	0.031–0.045
9	80.85–80.9	0.029–0.04
10	85.03–85.09	0.024–0.045
11	89.19–89.25	0.032–0.043
12	93.31–93.37	0.027–0.046
13	97.4–97.48	0.035–0.049
14	101.41–101.48	0.029–0.053

Allele (continued)	Mean	Standard Dev.
15	105.32–105.4	0.03–0.05
16	109.24–109.31	0.035–0.046
17	113.18–113.24	0.032–0.048
18	117.14–117.21	0.025–0.044
D12S391		
14	229.97–230.12	0.042–0.056
15	233.9–234.05	0.035–0.057
16	237.87–238.01	0.036–0.063
17	241.75–241.91	0.046–0.051
18	245.68–245.83	0.042–0.056
19	249.59–249.73	0.038–0.062
19.3	252.59–252.75	0.034–0.054
20	253.54–253.68	0.034–0.058
21	257.45–257.58	0.04–0.052
22	261.3–261.47	0.037–0.05
23	265.27–265.43	0.039–0.063
24	269.29–269.44	0.039–0.05
25	273.28–273.42	0.034–0.059
26	277.22–277.38	0.045–0.053
27	281.31–281.45	0.049–0.066
D16S539		
5	228.68–228.78	0.04–0.05
8	240.7–240.83	0.04–0.052
9	244.72–244.83	0.05–0.055
10	248.73–248.84	0.048–0.059
11	252.73–252.87	0.046–0.063
12	256.78–256.89	0.04–0.055
13	260.8–260.93	0.046–0.061
14	264.82–264.95	0.037–0.056
15	268.86–268.99	0.04–0.054
D18S51		
7	262.24–262.34	0.04–0.068
9	270.36–270.49	0.046–0.062
10	274.43–274.56	0.04–0.063
10.2	276.45–276.56	0.05–0.065
11	278.5–278.62	0.048–0.069
12	282.56–282.71	0.059–0.075
13	286.64–286.78	0.038–0.073

Allele <i>(continued)</i>	Mean	Standard Dev.
13.2	288.67–288.79	0.057–0.073
14	290.76–290.88	0.046–0.069
14.2	292.75–292.9	0.061–0.075
15	294.83–294.98	0.055–0.082
16	298.95–299.08	0.056–0.077
17	303.1–303.23	0.053–0.088
18	307.24–307.39	0.06–0.079
19	311.39–311.52	0.066–0.086
20	315.53–315.65	0.056–0.087
21	319.66–319.8	0.062–0.077
22	323.86–323.99	0.063–0.082
23	327.92–328.04	0.061–0.076
24	332.03–332.16	0.058–0.079
25	336.16–336.27	0.056–0.085
26	340.27–340.39	0.072–0.079
27	344.39–344.5	0.057–0.077
D19S433		
9	127.1–127.2	0.027–0.05
10	130.98–131.06	0.033–0.05
11	134.88–134.96	0.032–0.052
12	138.83–138.9	0.023–0.045
12.2	140.96–141.03	0.027–0.043
13	143.04–143.13	0.03–0.049
13.2	145.21–145.28	0.029–0.041
14	147.28–147.37	0.026–0.051
14.2	149.44–149.52	0.03–0.047
15	151.48–151.57	0.028–0.04
15.2	153.54–153.63	0.036–0.047
16	155.53–155.64	0.032–0.048
16.2	157.58–157.68	0.039–0.051
17	159.55–159.67	0.04–0.056
17.2	161.54–161.66	0.039–0.047
D1S1656		
9	175.77–175.96	0.046–0.09
10	179.67–179.85	0.047–0.08
11	183.53–183.71	0.054–0.08
12	187.39–187.57	0.056–0.078
13	191.24–191.4	0.04–0.069

Allele <i>(continued)</i>	Mean	Standard Dev.
14	195.13–195.29	0.04–0.062
14.3	198.01–198.16	0.04–0.061
15	198.87–199.02	0.045–0.07
15.3	201.86–202.01	0.046–0.062
16	202.72–202.88	0.034–0.06
16.3	205.76–205.91	0.05–0.06
17	206.63–206.79	0.044–0.063
17.3	209.66–209.82	0.048–0.064
18.3	213.56–213.73	0.047–0.061
19.3	217.48–217.64	0.044–0.073
20.3	221.4–221.56	0.042–0.064
D21S11		
24	185.02–185.14	0.042–0.069
24.2	187.02–187.12	0.042–0.066
25	188.98–189.08	0.045–0.066
26	192.92–193.01	0.037–0.056
27	196.87–196.96	0.036–0.052
28	200.75–200.83	0.032–0.052
28.2	202.71–202.81	0.036–0.053
29	204.69–204.78	0.035–0.045
29.2	206.73–206.82	0.039–0.049
30	208.71–208.8	0.04–0.048
30.2	210.68–210.78	0.039–0.049
31	212.7–212.79	0.034–0.05
31.2	214.67–214.76	0.037–0.049
32	216.69–216.78	0.04–0.046
32.2	218.65–218.75	0.04–0.048
33	220.67–220.76	0.031–0.056
33.2	222.62–222.72	0.036–0.059
34	224.74–224.85	0.039–0.048
34.2	226.65–226.74	0.042–0.055
35	228.7–228.8	0.042–0.057
35.2	230.64–230.73	0.041–0.052
36	232.62–232.74	0.042–0.046
37	236.68–236.78	0.037–0.052
38	240.61–240.71	0.04–0.052
D22S1045		
8	79.99–80.08	0.034–0.058

Allele (continued)	Mean	Standard Dev.
9	83.08–83.18	0.037–0.058
10	86.14–86.25	0.041–0.063
11	89.2–89.33	0.043–0.065
12	92.26–92.37	0.047–0.067
13	95.29–95.43	0.045–0.069
14	98.33–98.46	0.044–0.07
15	101.3–101.42	0.054–0.066
16	104.2–104.32	0.043–0.064
17	107.13–107.24	0.054–0.068
18	110.06–110.17	0.052–0.069
19	113.02–113.13	0.046–0.059
D2S1338		
15	289.59–289.7	0.038–0.054
16	293.68–293.78	0.036–0.053
17	297.73–297.82	0.042–0.055
18	301.81–301.9	0.046–0.064
19	305.9–306	0.043–0.06
20	310.03–310.11	0.047–0.057
21	314.12–314.22	0.042–0.062
22	318.22–318.31	0.034–0.06
23	322.3–322.39	0.044–0.062
24	326.38–326.46	0.044–0.058
25	330.46–330.55	0.038–0.06
26	334.54–334.61	0.047–0.056
27	338.63–338.71	0.046–0.057
28	342.95–343.05	0.037–0.061
D2S441		
9	80.29–80.38	0.028–0.054
10	84.41–84.5	0.033–0.057
11	88.51–88.62	0.03–0.062
12	91.66–91.76	0.046–0.067
13	92.58–92.69	0.045–0.062
14	96.49–96.62	0.036–0.069
15	100.51–100.63	0.045–0.067
16	104.4–104.53	0.038–0.056
D3S1358		
12	134.88–134.99	0.043–0.064
13	138.98–139.12	0.042–0.064

Allele <i>(continued)</i>	Mean	Standard Dev.
14	143.23–143.38	0.035–0.051
15	147.4–147.55	0.047–0.057
16	151.78–151.91	0.042–0.063
17	155.99–156.16	0.051–0.074
18	160.04–160.21	0.063–0.075
19	163.88–164.06	0.064–0.079
D8S1179		
8	123.54–123.64	0.04–0.068
9	127.61–127.71	0.039–0.071
10	131.7–131.81	0.045–0.064
11	135.86–135.95	0.045–0.07
12	140.1–140.2	0.046–0.056
13	144.67–144.77	0.043–0.058
14	149.13–149.23	0.046–0.07
15	153.45–153.57	0.053–0.069
16	157.7–157.85	0.054–0.073
17	161.86–162	0.056–0.086
18	165.94–166.07	0.054–0.085
19	169.99–170.15	0.054–0.096
FGA		
17	233.22–233.34	0.032–0.044
18	237.01–237.13	0.036–0.047
19	240.8–240.93	0.04–0.053
20	244.61–244.74	0.04–0.049
21	248.43–248.55	0.038–0.045
22	252.24–252.35	0.034–0.057
23	256.09–256.19	0.028–0.047
24	259.9–260.01	0.041–0.051
25	263.74–263.86	0.035–0.049
26	267.59–267.7	0.035–0.055
26.2	269.46–269.57	0.029–0.045
27	271.5–271.62	0.037–0.054
28	275.35–275.47	0.036–0.046
29	279.2–279.31	0.03–0.051
30	282.97–283.07	0.036–0.053
30.2	285.24–285.36	0.035–0.053
31.2	289.12–289.23	0.028–0.065
32.2	292.98–293.09	0.042–0.054

Allele (continued)	Mean	Standard Dev.
33.2	296.88–296.97	0.044–0.051
42.2	332.47–332.54	0.03–0.062
43.2	336.37–336.45	0.031–0.055
44.2	340.44–340.51	0.034–0.055
45.2	344.38–344.45	0.033–0.055
46.2	348.02–348.09	0.036–0.044
47.2	351.82–351.85	0.036–0.05
48.2	355.71–355.78	0.031–0.052
50.2	363.15–363.2	0.032–0.045
51.2	366.92–366.96	0.034–0.052
TH01		
4	181.53–181.67	0.044–0.065
5	185.44–185.58	0.039–0.063
6	189.32–189.46	0.037–0.066
7	193.2–193.33	0.028–0.058
8	197.05–197.17	0.032–0.053
9	200.93–201.03	0.036–0.047
9.3	203.93–204.03	0.037–0.052
10	204.85–204.95	0.031–0.051
11	208.78–208.9	0.036–0.048
13.3	219.64–219.75	0.038–0.047
vWA		
11	153.74–153.84	0.041–0.065
12	157.91–158.03	0.038–0.067
13	162.04–162.18	0.039–0.069
14	166.28–166.42	0.054–0.078
15	170.22–170.34	0.047–0.081
16	0.043–0.087	0.043–0.087
17	178.23–178.36	0.039–0.074
18	182.17–182.29	0.044–0.074
19	186.13–186.25	0.044–0.075
20	190.06–190.2	0.039–0.067
21	193.95–194.08	0.039–0.058
22	197.85–197.97	0.04–0.057
23	201.69–201.79	0.032–0.051
24	206.01–206.12	0.032–0.063

Extra peaks in the electropherogram

Causes of extra peaks

Peaks other than the target alleles may be detected on the electropherogram. Causes for the appearance of extra peaks include stutter products, incomplete 3' A nucleotide addition (at the n-1 position), dye artifacts, and mixed DNA samples (see DAB Standard 8.1.2.2).

Stutter products

Stutter is a well-characterized PCR artifact that refers to the appearance of a minor peak one repeat unit smaller (or less frequently, one repeat larger) than the major STR product (Butler, 2005; Mulero *et al.*, 2006). 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. Peak heights were measured for amplified samples (n = 1023) at the loci used in the NGM™ Kit. All data were generated on the Applied Biosystems™ 3130xl Genetic Analyzer.

Due to the enhanced buffer system developed for the NGM™ Kit, plus stutter peaks may be encountered at other loci more frequently than with older AmpFSTR™ chemistries (for example, the SGM Plus™ Kit). Laboratories should evaluate the occurrence of plus stutter peaks and develop appropriate interpretation guidelines as part of an internal validation study.

Some conclusions from these measurements and observations are:

- For each NGM™ Kit locus, the percent stutter generally increases with allele length, as shown in Figure 9 to Figure 13 on page 73 to 75.
- Smaller alleles display a lower level of stutter relative to the longer alleles within each locus.
- Each allele within a locus displays a percent stutter that is consistent.
- Stutter filter sets in GeneMapper™ ID Software and GeneMapper™ ID-X Software, calculated as the mean stutter for the locus plus three standard deviations (n = 1023), are shown in Table 5 on page 75. Peaks in the stutter position that are above the stutter filter percentage specified in the software are not filtered. Peaks in the stutter position that have not been filtered and remain labeled can be further evaluated. For evaluation of mixed samples, see Figure 24 on page 90.
- The measurement of percent stutter for allele peaks that are off-scale may be unusually high due to artificial truncation of the main allele peak.

Figure 9 Stutter percentages for D10S1248, D12S391 and D16S539 loci (Blue and red colors indicate loci labeled with FAM™ and PET™ dyes respectively)

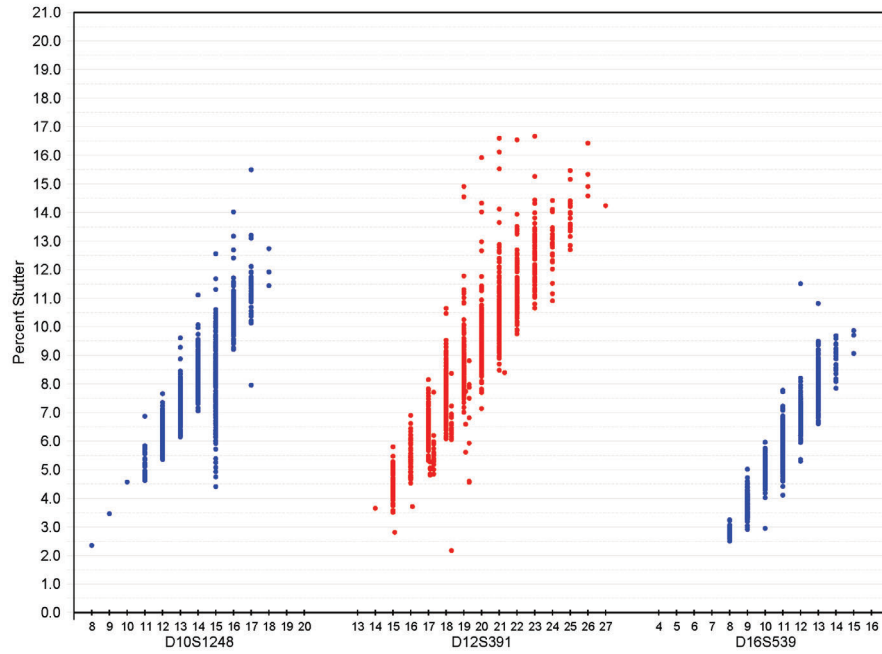


Figure 10 Stutter percentages for the D18S51, D19S433 and D1S1656 loci (Green, black, and red colors indicate loci labeled with VIC™, NED™, and PET™ dyes, respectively)

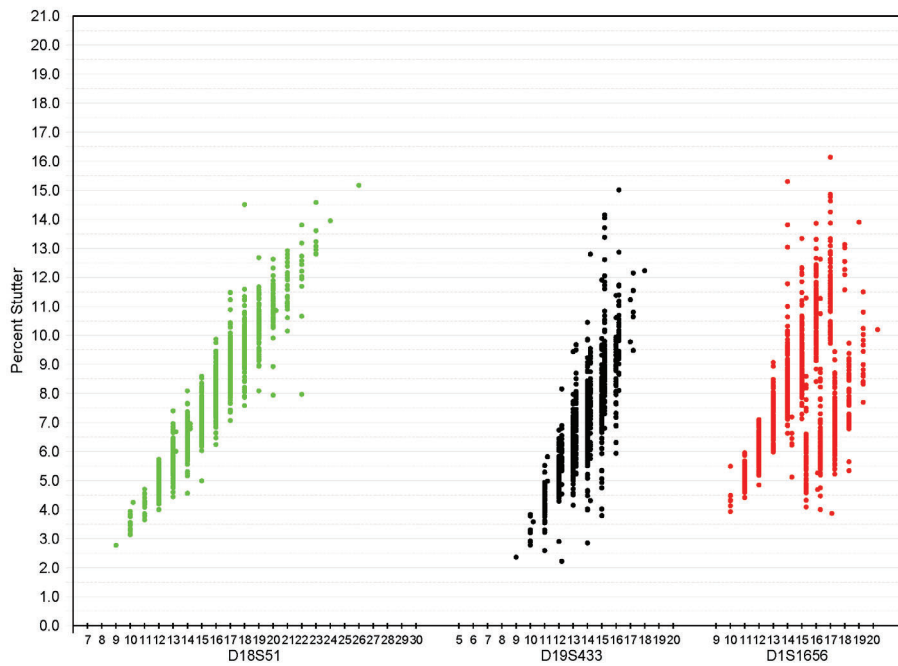


Figure 11 Stutter percentages for the D21S11, D22S1045 and D2S1338 loci (Green, black/gray, and blue colors indicate loci labeled with VIC™, NED™ and FAM™ dyes, respectively. Black and gray data points associated with the D22S1045 locus indicate minus- and plus-stutter, respectively)

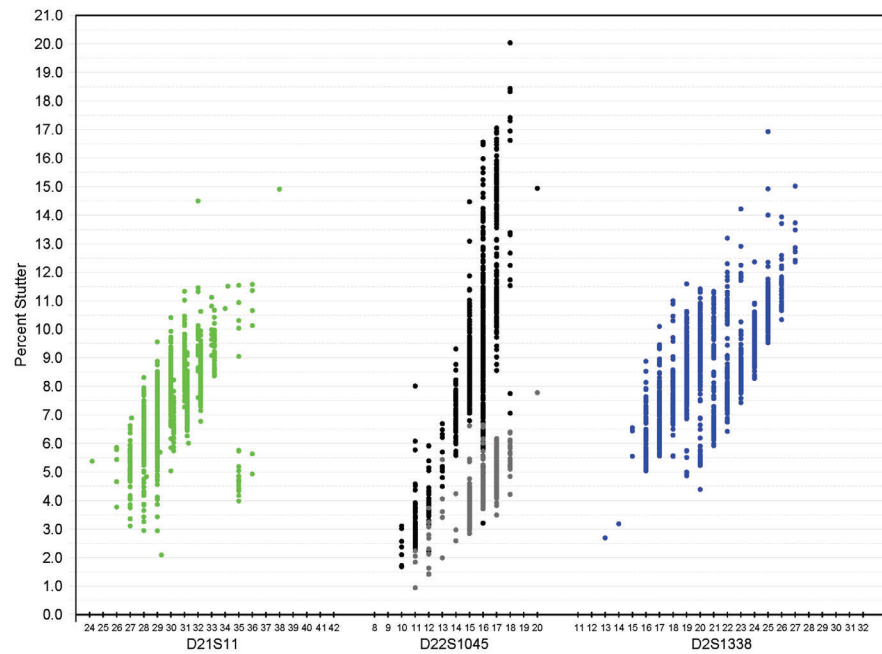


Figure 12 Stutter percentages for the FGA and TH01 loci (Black data points indicate loci labeled with NED™ dye)

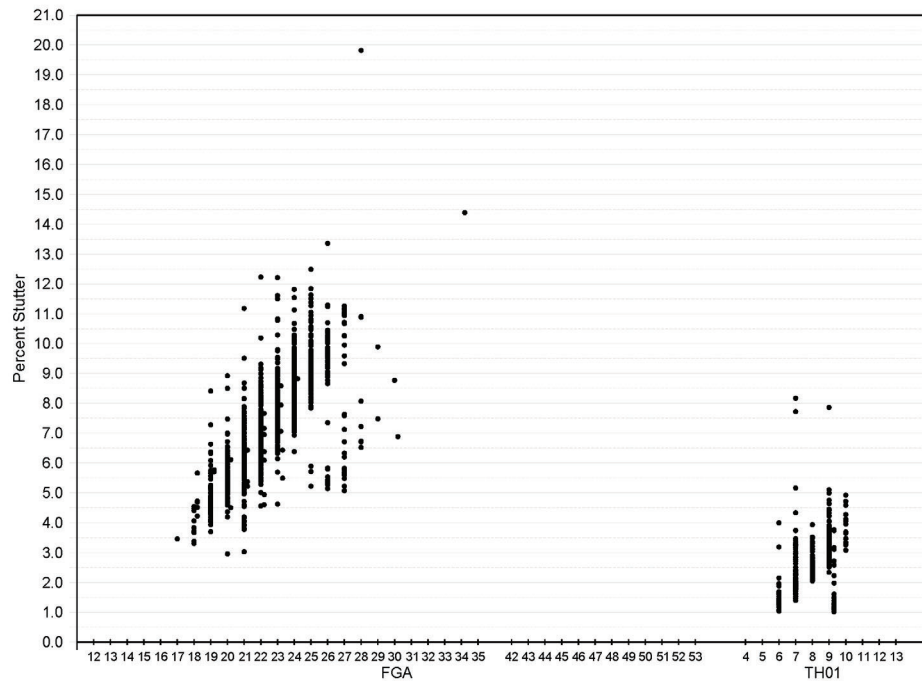


Figure 13 Stutter percentages for the D2S441, D3S1358, D8S1179, and vWA loci (Red, green, and blue colors indicate loci labeled with and PET™, VIC™, and FAM™ dyes, respectively)

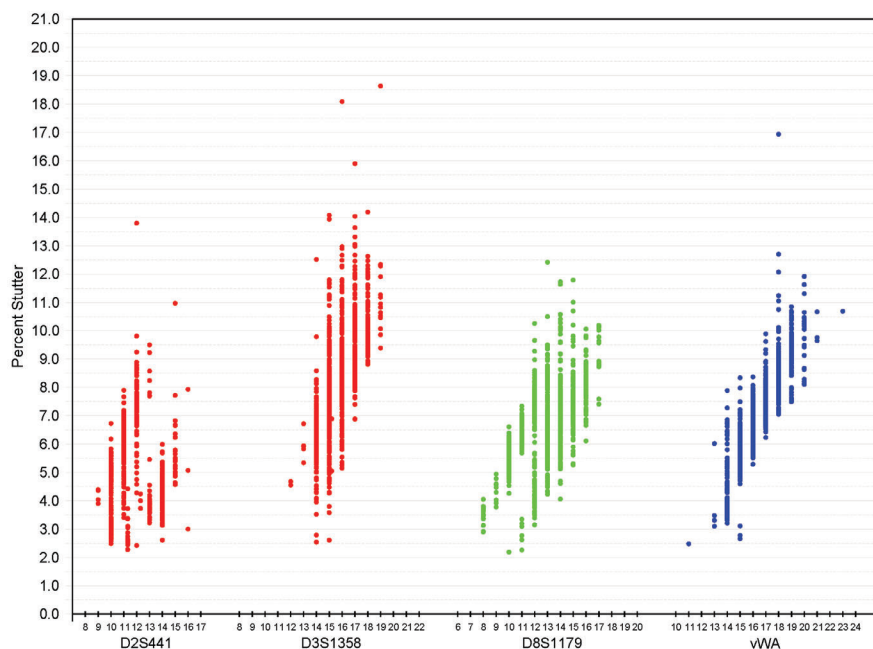


Table 5 Marker-specific stutter filter percentages for NGM™ Kit loci

Locus	% Stutter
D10S1248	12.39
vWA	11.83
D16S539	10.12
D2S1338	12.83
D8S1179	10.31
D21S11	10.87
D18S51	14.08
D22S1045 (-3nt)	18.05
D22S1045 (+3 nt)	7.63
D19S433	11.20
TH01	4.27
FGA	12.10
D2S441	9.45
D3S1358	13.07
D1S1656 (-4nt)	14.46
D1S1656 (-2nt) [†]	4.70
D12S391	15.27

[†] The - 2nt stutter filter is not included in GeneMapper™ ID Software v3.2.1NGM_panel_v2 due to functional limitations of the software.

IMPORTANT! The values shown are the values we determined during developmental validation studies. We recommend that laboratories perform internal validation studies to determine the appropriate values to use.

Addition of 3' A nucleotide

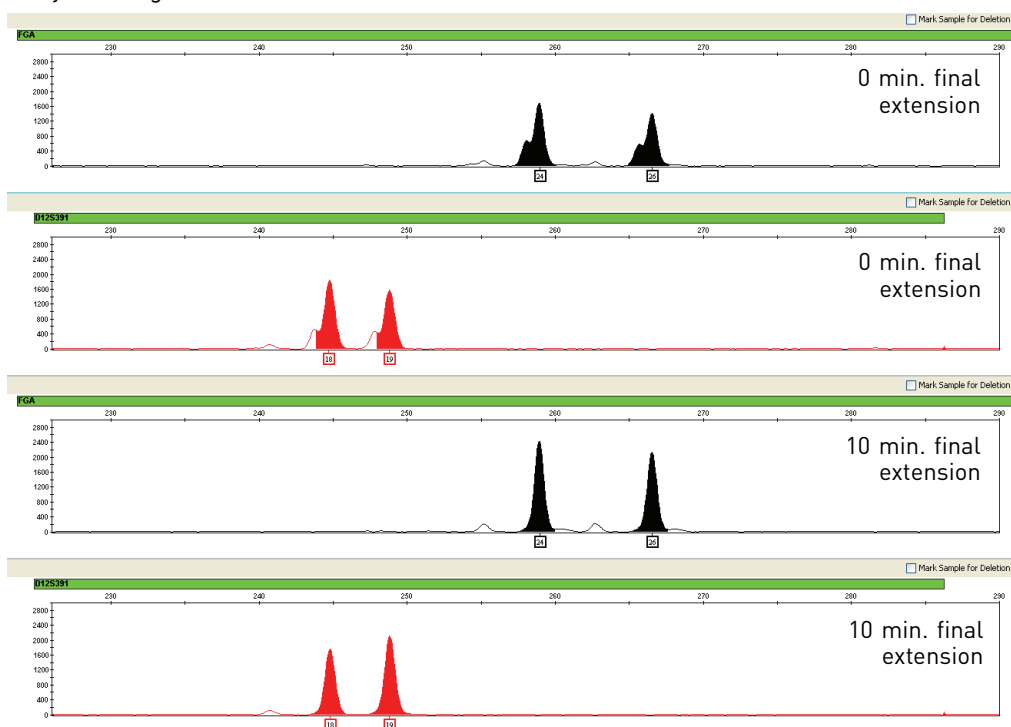
Many DNA polymerases can catalyze the addition of a single nucleotide (predominantly adenosine) to the 3' ends of double-stranded PCR products (Clark, 1988; Magnuson *et al.*, 1996). This non-template addition results in a PCR product that is one nucleotide longer than the actual target sequence. The 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 NGM™ Kit includes two main design features that promote maximum +A addition:

- The primer sequences have been optimized to encourage +A addition.
- The new, highly robust PCR chemistry allows complete +A addition with a short final incubation at 60 °C for 10 minutes.

This final extension step gives the DNA polymerase additional time to complete +A addition to all double-stranded PCR products. See Figure 14 on page 76 for examples of incomplete and normal +A addition. Final extension incubation for longer than the recommended 10 minutes may result in double +A addition, in which *two* non-template adenosine residues are added to the PCR product. Double +A addition can cause "shoulders" on the right side of main allele peaks, and is therefore to be avoided.

Figure 14 Omitting the final extension step results in shoulders on main allele peaks due to incomplete A nucleotide addition. Data are from an Applied Biosystems™ 3130xl Genetic Analyzer using the NGM™ Kit



Due to improved PCR buffer chemistry, the lack of +A addition is generally less an issue with the NGM™ Kit than with earlier generation kits. However, “shouldering” of allele peaks may still be observed if the amount of input DNA is greater than that recommended by the NGM™ Kit protocol. Amplification of excess input DNA may also result in offscale data.

Artifacts

Artifacts and anomalies are seen in all molecular biological systems. Artifacts are typically reproducible while anomalies are non-reproducible, intermittent occurrences that are not observed consistently in a system (for example, spikes and baseline noise). Due to improvements in PCR primer manufacturing processes, the incidence of artifacts has been greatly reduced in the NGM™ Kit. NGM™ Kit electropherograms are essentially free of reproducible dye artifacts within the Kit's read region of 68-407 nt. Figure 15 on page 78 shows the very low baseline level fluorescence of a typical negative control PCR using the NGM™ Kit.

Most STR loci produce minus-stutter peaks as a by-product of PCR amplification. A process of “slippage” has been proposed as a molecular mechanism for stutter, where the Taq DNA polymerase enzyme “slips” on the template DNA during replication and produces a minority PCR product that is shorter than the template strand, usually by one repeat unit. The stutter process may also occur in the opposite direction to produce amplicon DNA that is usually one repeat unit longer than the template strand, termed plus-stutter. While plus-stutter is normally much less significant than minus-stutter in STR loci with tetranucleotide repeats, the incidence of plus-stutter may be more significant in trinucleotide repeat-containing loci. The D22S1045 locus in the NGM™ Kit is a trinucleotide repeat locus, and shows an elevated level of plus-stutter. For example, Figure 17 on page 79 is an electropherogram of the D22S1045 locus showing plus stutter. GeneMapper™ ID Software and GeneMapper™ ID-X Software analysis parameter files supplied for use with the NGM™ Kit contain a plus-stutter filter for the D22S1045 locus to prevent these peaks from being called in normal profiles.

Plus stutter may also be visible at tetranucleotide repeat loci in next generation NGM™ Kits due to improvements in buffer formulation over previous kits. Laboratories should evaluate the occurrence of plus stutter across the profile as part of internal validation studies.

Figure 16 on page 78 shows an example of a non-standard (minus 2-nt) stutter that may be observed in certain STR loci such as D1S1656 with more complex nucleotide sequences that include regions of dinucleotide TG repeats. Genotyping may result in the detection of these artifacts as off-ladder (OL) alleles.

It is important to consider possible noise and artifacts when interpreting data from the NGM™ Kit on the Applied Biosystems™ 3500/3500xL and 3130/3130xL, ABI PRISM™ 3100/3100-*Avant*, and ABI PRISM™ 310 Genetic Analyzers. Note that a high degree of magnification is used in the sample electropherograms shown in Figure 15 to Figure 17 on page 78 to 79.

Figure 15 Examples of fluorescence background in data produced on an Applied Biosystems™ 3130xl Genetic Analyzer, (Y-axis scale 0–100 RFU)

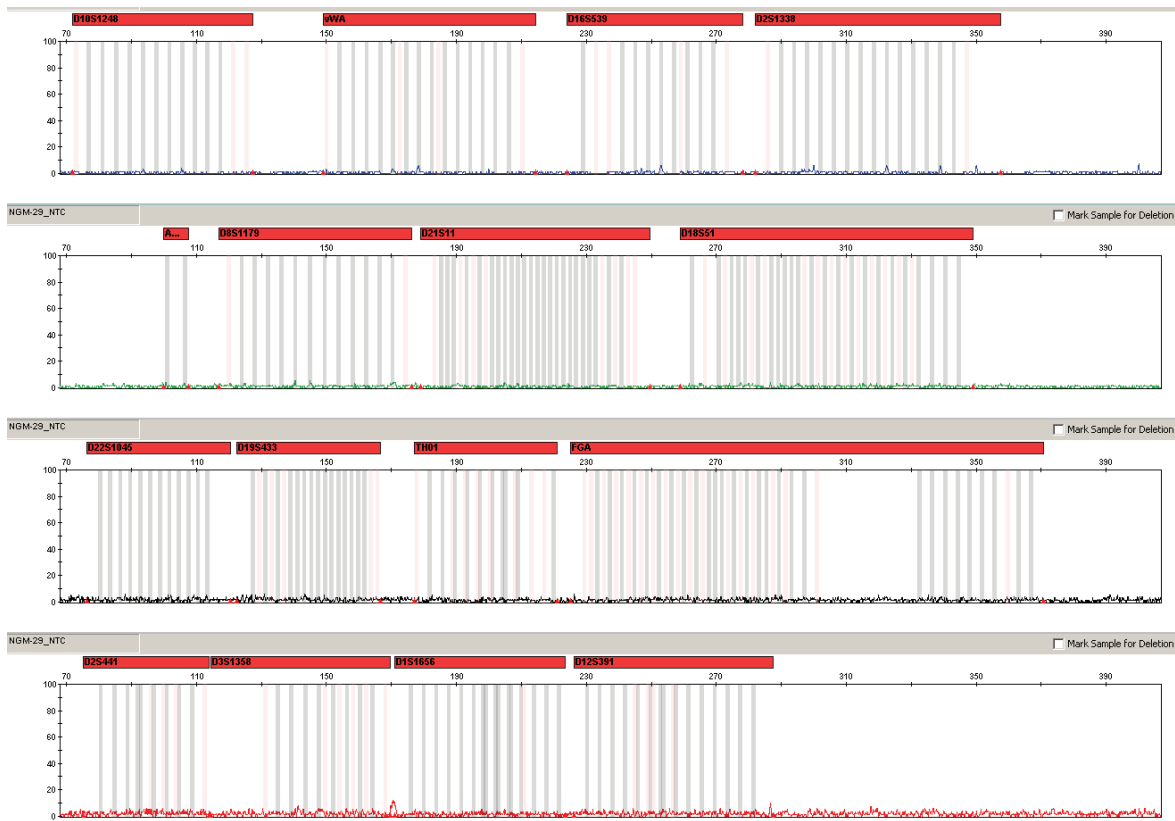


Figure 16 Example of a -2 nt reproducible artifact at the D1S1656 locus. Data produced on an Applied Biosystems™ 3130xl Genetic Analyzer

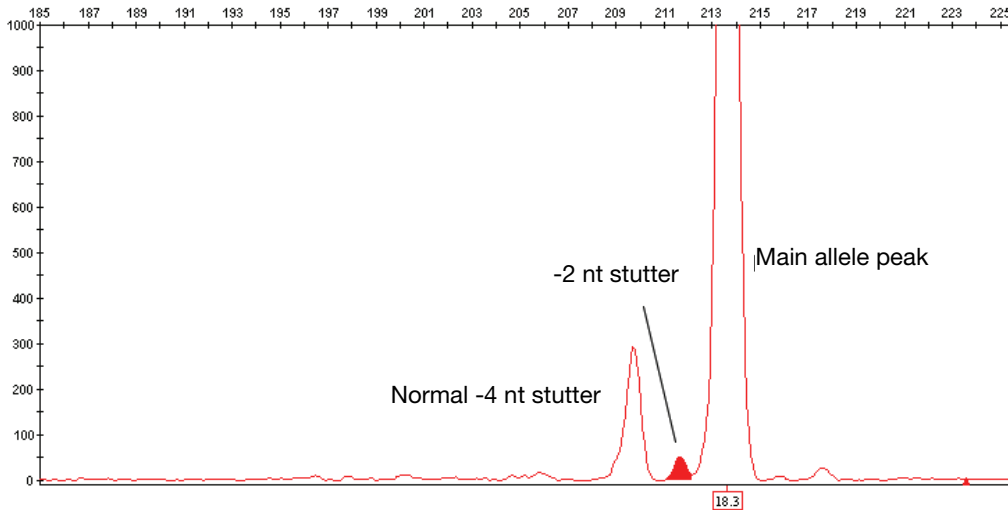
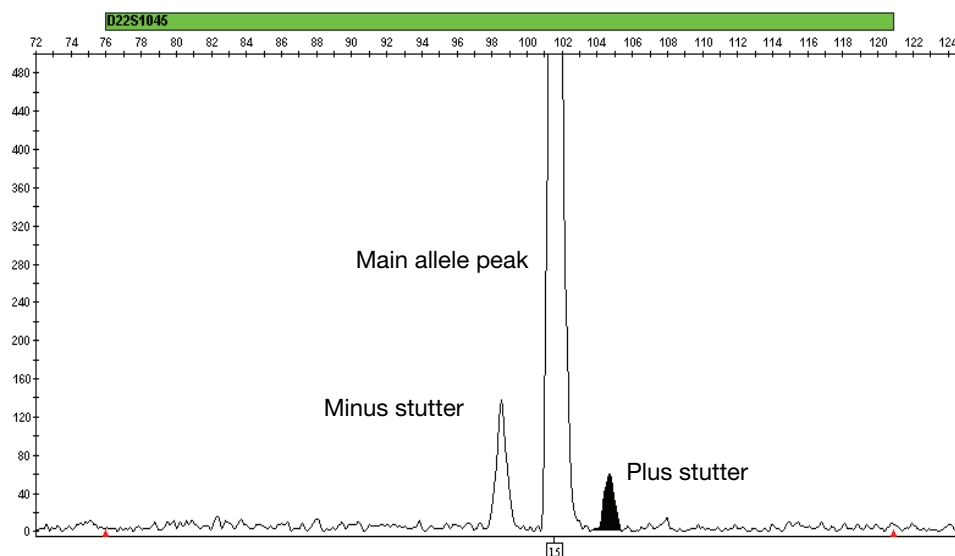


Figure 17 NGM™ Kit electropherogram showing plus stutter associated with the D22S1045 STR locus. Data produced on an Applied Biosystems™ 3130xl Genetic Analyzer



Characterization of loci

SWGAM guideline 2.1

“The basic characteristics of a genetic marker must be determined and documented.”
(SWGAM, July 2003)

This section describes basic characteristics of the 15 loci and the sex-determining marker, Amelogenin, which are amplified with the NGM™ Kit. Most of these loci have been extensively characterized by other laboratories.

Nature of the polymorphisms

The primers for the Amelogenin locus flank a 6-nucleotide deletion within intron 1 of the X homolog. Amplification results in 104-nt and 110-nt products from the X and Y chromosomes, respectively. (Sizes are the actual nucleotide size according to sequencing results, including 3' A nucleotide addition, and size may not correspond exactly to allele mobility observed on capillary electrophoresis platforms.) With the sole exception of D22S1045, a trinucleotide STR, the remaining NGM™ Kit loci are tetranucleotide short tandem repeat (STR) loci. The length differences among alleles of a particular locus result from differences in the number of repeat units.

All the alleles in the AmpF ϕ STR™ NGM™ Allelic Ladder, including microvariants, have been subjected to sequencing at Life Technologies. In addition, other groups in the scientific 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 NGM™ Kit loci, there is consensus on the repeat patterns and structure of the STRs.

Inheritance

The Centre d'Etude du Polymorphisme Humain (CEPH) has collected DNA from 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).

Mapping

The NGM™ Kit loci 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).

Genetic linkage

Two sets of STR loci in the NGM™ Kit are located on the same chromosomes. vWA and D12S391 are located approximately 6.3 million bp apart on the p arm of chromosome 12, while D2S1338 and D2S441 are located approximately 150 million bp apart on opposite arms of chromosome 2. Linkage disequilibrium analysis was conducted on the genotype results from 996 individuals of three ethnic groups (see "Population data" on page 91). STR locus genotype results from the population study were analyzed using the Linkage Disequilibrium module of GenePop software version 4.0.10 (Raymond and Rousset, 1995; Rousset, 2008). See Table 6 for results.

The relatively high probability values indicate that there is no statistically significant linkage disequilibrium found between the pairs of loci located on the same chromosome.

An independent analysis of the same set of population data by Budowle *et al.* (2010) concluded that there was no evidence of LD between or within any of the NGM™ Kit, and that the multiplication rule could therefore be used for all markers contained in the 15 STR markers contained in the kit for the purpose of calculating the rarity of a DNA profile. However, they cautioned that, while there was no evidence of LD at the population level, the independence of vWA and D12S391 could not be assumed for the purpose of kinship analysis, due to the proximity of the loci on Chromosome 12.

Table 6 GenePop LD Result (p value for pairwise analysis of loci)

Locus	Chromosome Map Position†	Chromosome Nuclear Coordinates† (million bp)	African-American (N = 350)	Caucasian (N = 350)	Hispanic (N = 293)
vWA	p13.31	5.9	0.86	0.29	0.27
D12S391	p13.2	12.2			
D2S441	p14	68	0.11	0.32	0.19
D2S1338	q35	218			

† STR locus mapping data was obtained from the NCBI Map Viewer http://www.ncbi.nlm.nih.gov/projects/mapview/map_search.cgi?taxid=9606 or the UCSC Genome Browser (<http://genome.ucsc.edu/>). GenePop LD analysis probability results (p values) greater than 0.05 were considered to indicate that linkage disequilibrium between the loci within the population tested was not statistically significant.

Species specificity

SWGAM Guideline 2.2

“For techniques designed to type human DNA, the potential to detect DNA from forensically relevant nonhuman species should be evaluated.” (SWGAM, July 2003)

The NGM™ Kit provides the required specificity for detecting human alleles.

Nonhuman studies

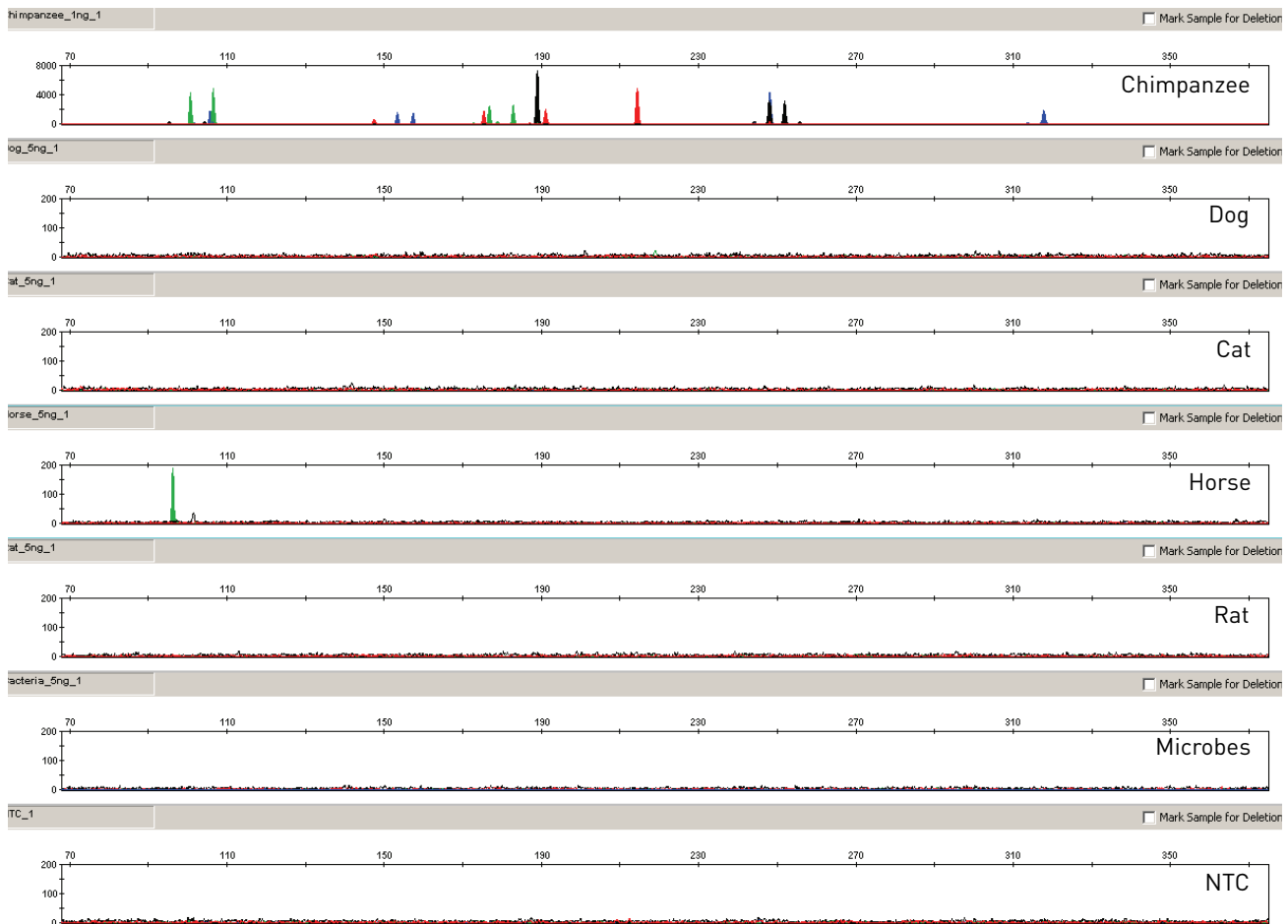
Nonhuman DNA may be present in forensic casework samples. The following species were tested (in the specified amounts) using standard PCR and capillary electrophoresis conditions for the NGM™ Kit.

- Primates: gorilla, chimpanzee, and macaque (1.0 ng each)
- Non-primates: mouse, dog, sheep, pig, rabbit, cat, horse, hamster, rat, chicken, and cow (5.0 ng each)
- Microorganisms: *Candida albicans*, *Staphylococcus aureus*, *Escherichia coli*, *Neisseria gonorrhoeae*, *Bacillus subtilis*, and *Lactobacillus rhamnosus* (equivalent to 105 copies)

Results were assessed for the presence of any amplified peaks that would indicate cross reactivity of the NGM™ Kit with any of these non-human species.

Figure 18 on page 82 shows example electropherogram results from the species specificity tests. The chimpanzee and gorilla DNA samples produced partial profiles within the 70 – 283 nucleotide region (gorilla data not shown). Macaque DNA produced a strong Amelogenin-X peak and two small out-of-marker-range peaks in PET (data not shown).

Figure 18 Representative electropherograms from a species specificity study including positive and non-template controls (NTC)



The microorganisms, cow, sheep, pig, dog, cat, chicken, hamster, mouse, rabbit, and rat did not yield detectable products. Of the non-primates, only horse DNA produced a 96-bp fragment near the amelogenin locus in the VIC™ dye.

Sensitivity

SWGDM guideline 2.3

“When appropriate, the range of DNA quantities able to produce reliable typing results should be determined.” (SWGDM, July 2003)

Importance of quantification

The recommended amount of input DNA for the NGM™ Kit is 1.0 ng, based on quantification using either the Quantifiler™ Human or Quantifiler™ Duo Quantification Kit and individual laboratories should determine the optimum input DNA amount according to the quantification method in use in the laboratory. If the sample contains degraded or inhibited DNA, amplification of a higher concentration

of DNA may be beneficial. In Figure 19 on page 84, the control DNA 007 was serially diluted from 2.0 ng to 0.062 ng. Full profiles (32 PCR products) were consistently obtained at 0.125 ng, but occasional partial profiles missing 1 – 3 alleles were observed at 0.062 ng.

Effect of DNA quantity on results

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

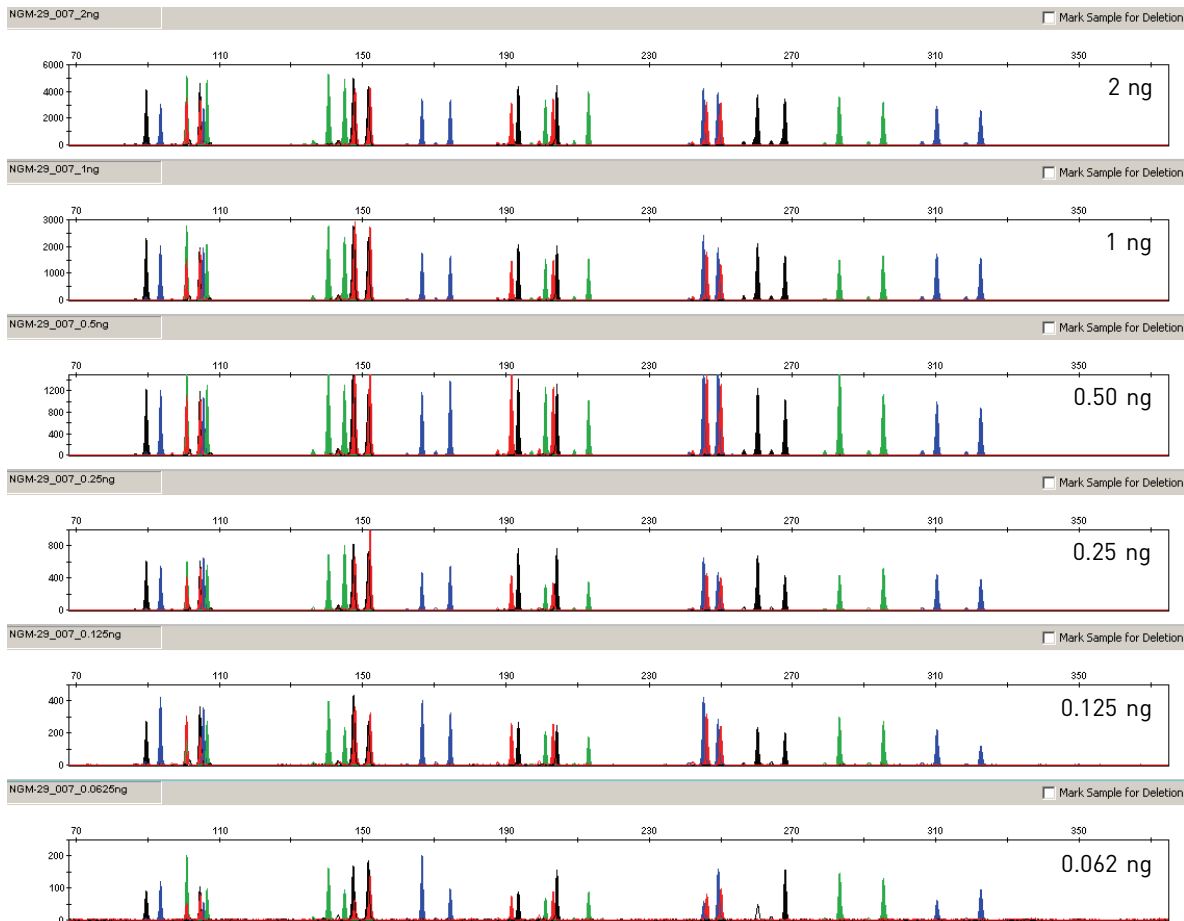
- Fluorescence intensity that exceeds the linear dynamic range for detection by the instrument (“off-scale” data).
- Off-scale data. Off-scale data is a problem because:
 - Quantification (peak height and area) for off-scale peaks is not accurate. For example, an allele peak that is off-scale can cause 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. This inaccuracy results in poor spectral separation (“pull-up”).
- Incomplete +A nucleotide addition.

To address these issues, reamplify the sample using less DNA.

When the total number of allele copies added to the PCR is extremely low, unbalanced amplification of the alleles may occur because of stochastic fluctuation.

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

Figure 19 Electropherograms for 29-cycle amplifications using 2 ng, 1 ng, 0.50 ng, 0.25 ng, 0.125 ng, and 0.062 ng of control DNA 007. Electrophoresis was performed on an Applied Biosystems™ 3130xl Genetic Analyzer. Note that the y-axis scale is magnified for the smaller input amounts of DNA



Stability

SWGDM guideline 2.4

“The ability to obtain results from DNA recovered from biological samples deposited on various substrates and subjected to various environmental and chemical insults has been extensively documented. In most instances, assessment of the effects of these factors on new forensic DNA procedures is not required. However, if substrates and/or environmental and/or chemical insults could potentially affect the analytical process, then the process should be evaluated using known samples to determine the effects of such factors.” (SWGDM, July 2003)

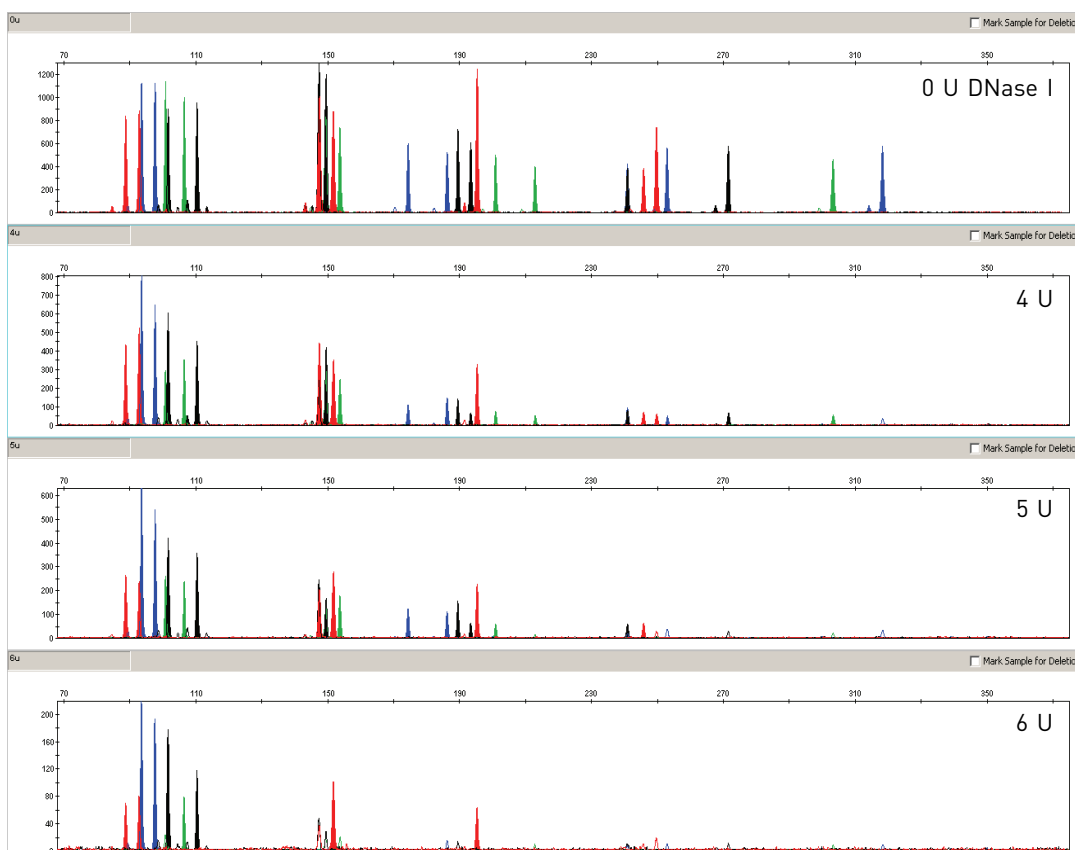
Degraded DNA

As the average size of degraded DNA approaches the size of the target sequence, the amount of PCR product generated is reduced because of 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 Raji DNA was sonicated and incubated with increasing doses of DNase I (0 to 6 Units) for 20 minutes (Bender *et al.*, 2004). The DNA was examined by agarose gel analysis to determine the average size of the DNA fragments at each time point.

Amplification of 1.0 ng of degraded DNA using the NGM™ Kit was performed. As the DNA became progressively degraded, the loci failed to amplify robustly in order of decreasing size. Preferential amplification was not observed.

Figure 20 Amplification of Raji DNA samples sonicated and incubated with increasing doses of DNase I. PCR amplification was done for 29 cycles, with electrophoresis being performed on an Applied Biosystems™ 3130xl. Panels 1, 2, 3, and 4 correspond to 0, 4, 5, and 6 units of DNase I. Note that the y-axis scale is magnified for more degraded samples, which generate lower peak heights



The larger loci contained in the NGM™ Kit, which fail to amplify in significantly degraded samples may be recovered by using the AmpF!STR™ MiniFler™ Kit. The amplicon size for the larger loci has been reduced to facilitate performance on degraded samples. For more information please refer to the *AmpF!STR™ MiniFler™ PCR Amplification Kit User's Guide* (Part no. 4374618).

Effect of inhibitors — hematin

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

To examine the effects of hematin on the performance of the NGM™ Kit, 1.0 ng of control DNA 007 was amplified in the presence of increasing concentrations of hematin for 29 cycles of amplification (Figure 21 on page 86). The concentrations of hematin used were 0 μM, 50 μM, 100 μM, 150 μM, and 200 μM (see Table 7 on page 86).

Figure 21 Electropherograms for the AmpF ϕ STR™ NGM™ and SGM Plus™ Kits show the improved performance of the NGM™ Kit in the presence of hematin compared with previous AmpF ϕ STR™ kits. In order from top to bottom, the panels show profiles for: NGM™ Kit uninhibited control, NGM™ Kit with 200 μM hematin, SGM Plus™ Kit uninhibited control and SGM Plus™ Kit with 200 μM hematin. The SGM Plus™ and NGM™ Kits were amplified for 28 and 29 cycles, respectively

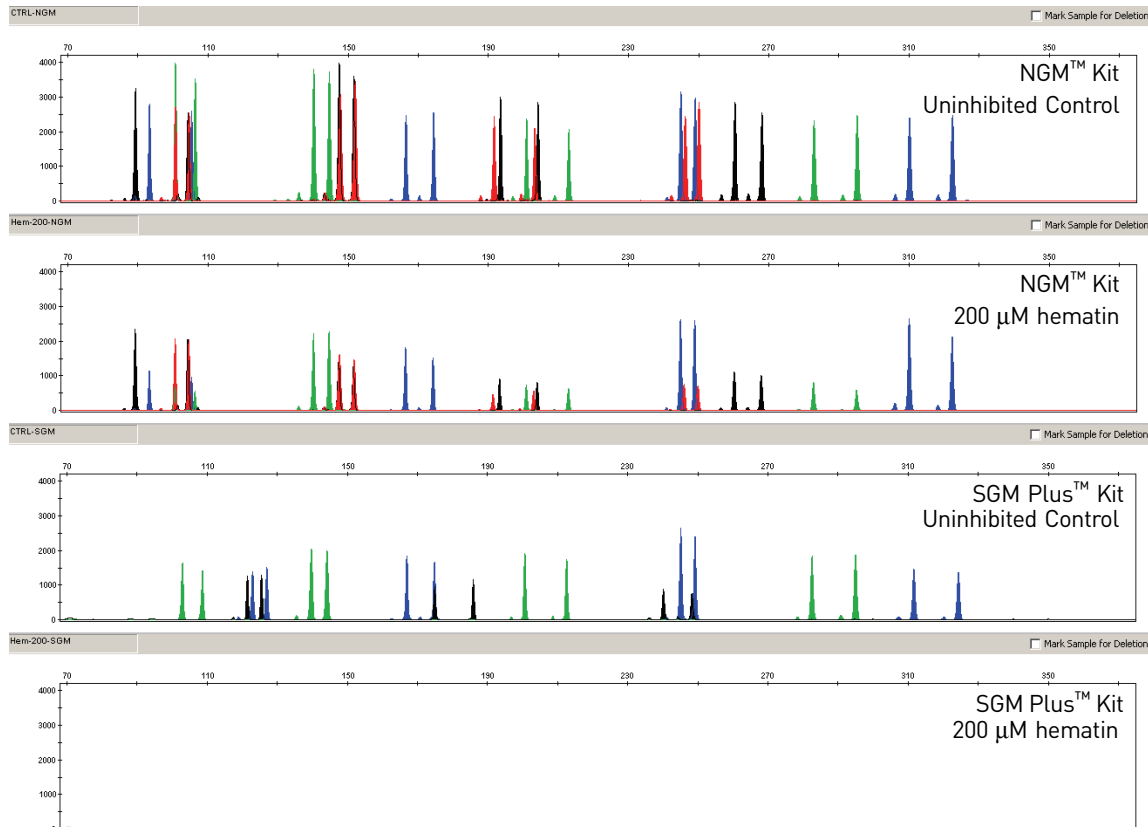


Table 7 NGM™ Kit performance in simulated hematin inhibition (n = 3)

Hematin Concentration	Number of Alleles Detected [†]
0 μM	32, 32, 32
50 μM	32, 32, 32
100 μM	32, 32, 32
150 μM	32, 32, 32
200 μM	32, 32, 32

[†] Only those peaks >50 RFU were counted. A complete profile with control DNA 007 yields 32 peaks using the NGM™ Kit on an Applied Biosystems™ 3130xl instrument.

Effect of inhibitors — humic acid

Traces of humic acid may inhibit the PCR amplification of DNA evidence collected from soil. Amplification of 1 ng of control DNA 007 in the presence of increasing amounts of humic acid was performed using the NGM™ Kit for 29 cycles of amplification (see Figure 22). The concentrations of humic acid tested were 0, 20, 40, 60, and 80 ng/μL (see Table 8 on page 87).

Figure 22 Electropherograms for the AmpF λ STR™ NGM™ and SGM Plus™ Kits show improved performance of the NGM™ Kit in the presence of humic acid compared to previous AmpF λ STR™ kits. In order from top to bottom, the panels show profiles for: NGM™ Kit uninhibited control, NGM™ Kit with 80 ng/μL humic acid, SGM Plus™ Kit uninhibited control, and SGM Plus™ Kit with 80 ng/μL humic acid. The SGM Plus™ and NGM™ Kits were amplified for 28 and 29 cycles, respectively

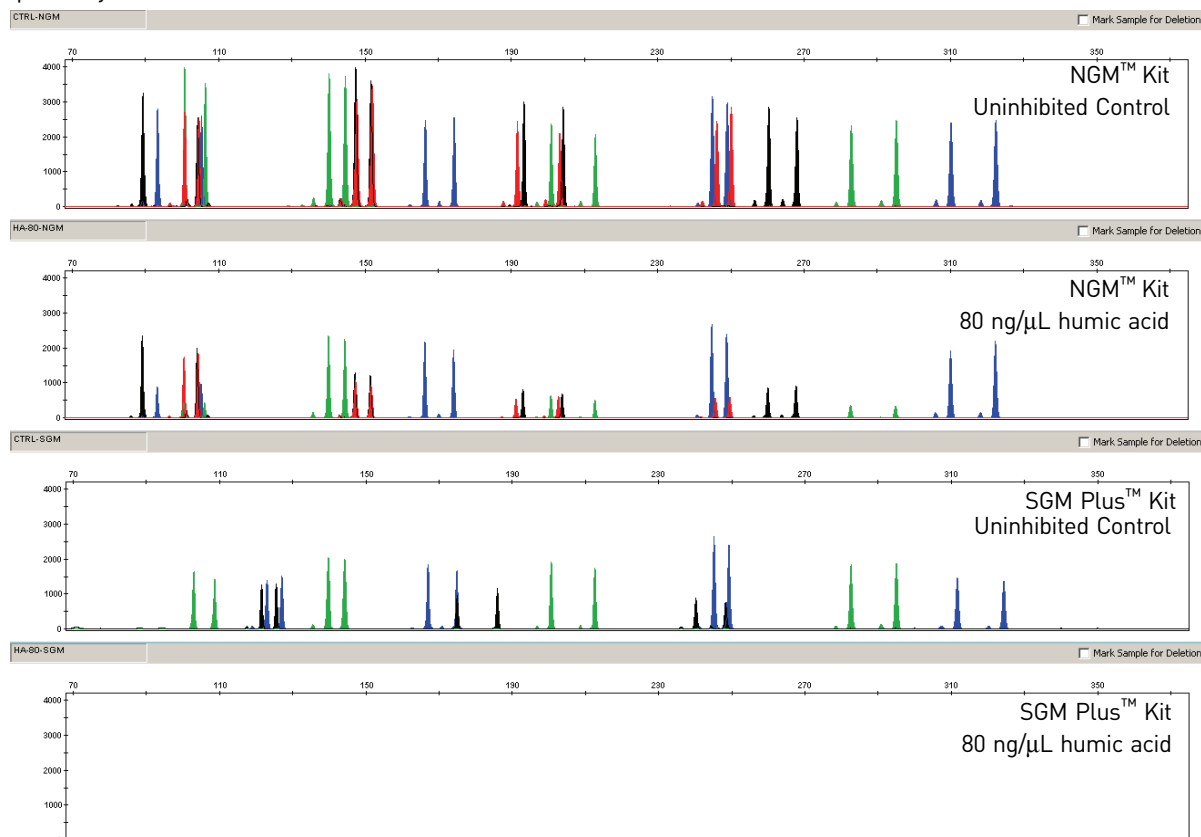


Table 8 NGM™ Kit performance in simulated model of humic acid inhibition (n = 3)

Humic acid concentration	Number of alleles detected [†]
0 ng/μL	32, 32, 32
20 ng/μL	32, 32, 32
40 ng/μL	32, 32, 32
60 ng/μL	32, 32, 32
80 ng/μL	32, 32, 32

[†] Only those peaks >50 RFU were counted. A complete profile with control DNA 007 yields 32 peaks using the NGM™ Kit on an Applied Biosystems™ 3130xl instrument.

Mixture studies

SWGDM guideline 2.8

“The ability to obtain reliable results from mixed source samples should be determined.” (SWGDM, July 2003)

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 determine 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.

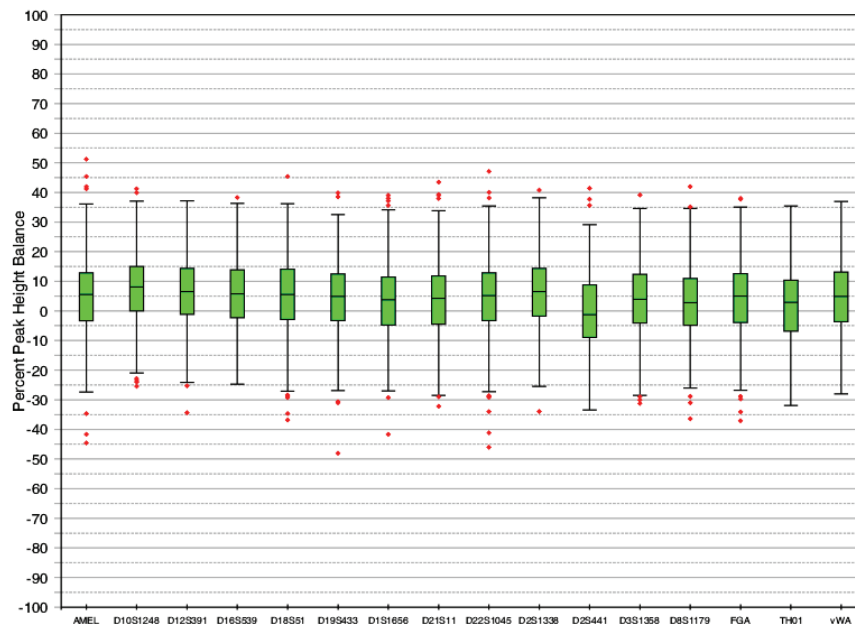
Mixture studies

Evidence samples that contain body fluids and/or tissues originating from more than one individual are an integral component of forensic casework. Therefore, it is essential to ensure that the DNA typing system is able to detect DNA mixtures. Mixed samples can be distinguished from single-source samples by:

- 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 typically observed in a single-source sample
- 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, median, minimum, and maximum peak height ratios observed for alleles in the NGM™ Kit loci in unmixed population database samples are shown in Figure 23 below.

Figure 23 Heterozygote ratios for 1 ng of input DNA amplified for 29 cycles with the NGM™ Kit. The distribution of intra-locus peak height ratios are expressed as plus and minus percent, by locus. Green boxes show the middle 50% or interquartile range (IQR). Box halves below and above median show the second and third quartile, respectively. “Whiskers” indicate 1.5 IQR from the upper and lower margins of the IQR. Red diamonds are outlier data points more than 1.5 IQR from the median



If an unusually low peak height ratio is observed 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. Possible causes of imbalance at a locus are:

- Degraded DNA
- Presence of inhibitors
- Extremely low amounts of input DNA
- A mutation in one of the primer binding sites
- Presence of an allele containing a rare sequence that does not amplify as efficiently as the other allele

Resolution of genotypes in mixed samples

A sample containing DNA from two sources can comprise (at a single locus) any of the seven genotype combinations (see 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 or not 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 Applied Biosystems™ instruments provides additional valuable data to aid in resolving mixed genotypes.

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 (0:1, 1:1, 3:1, 7:1, 15:1, 1:0). The total amount of genomic input DNA mixed at each ratio was 1.0 ng. The samples were amplified in a GeneAmp™ PCR System 9700, then electrophoresed and detected using an Applied Biosystems™ 3130xl Genetic Analyzer.

The results of the mixed DNA samples are shown in Figure 24 on page 90 where samples A and B were mixed according to the ratios indicated. The minor component allele calls at non-overlapping loci are highlighted. Detection of full profiles for the minor contributor was possible at ratios of 3:1 (0.750:0.250 ng) and 7:1 (0.875:0.125 ng). Generally, 15:1 ratios resulted in partial profiles for the minor component. The profiles of these samples are described in Table 9 on page 90.

Figure 24 Amplification of DNA mixtures at various ratios. Panels show electropherograms for (top to bottom): Major contributor only, 1:1 mixture (maj:min), 3:1 mixture, 7:1 mixture, 10:1 mixture and 15:1 mixture. The experiment was performed with both 29- and 30 cycle amplification; electropherograms shown are from the 29-cycle amplification

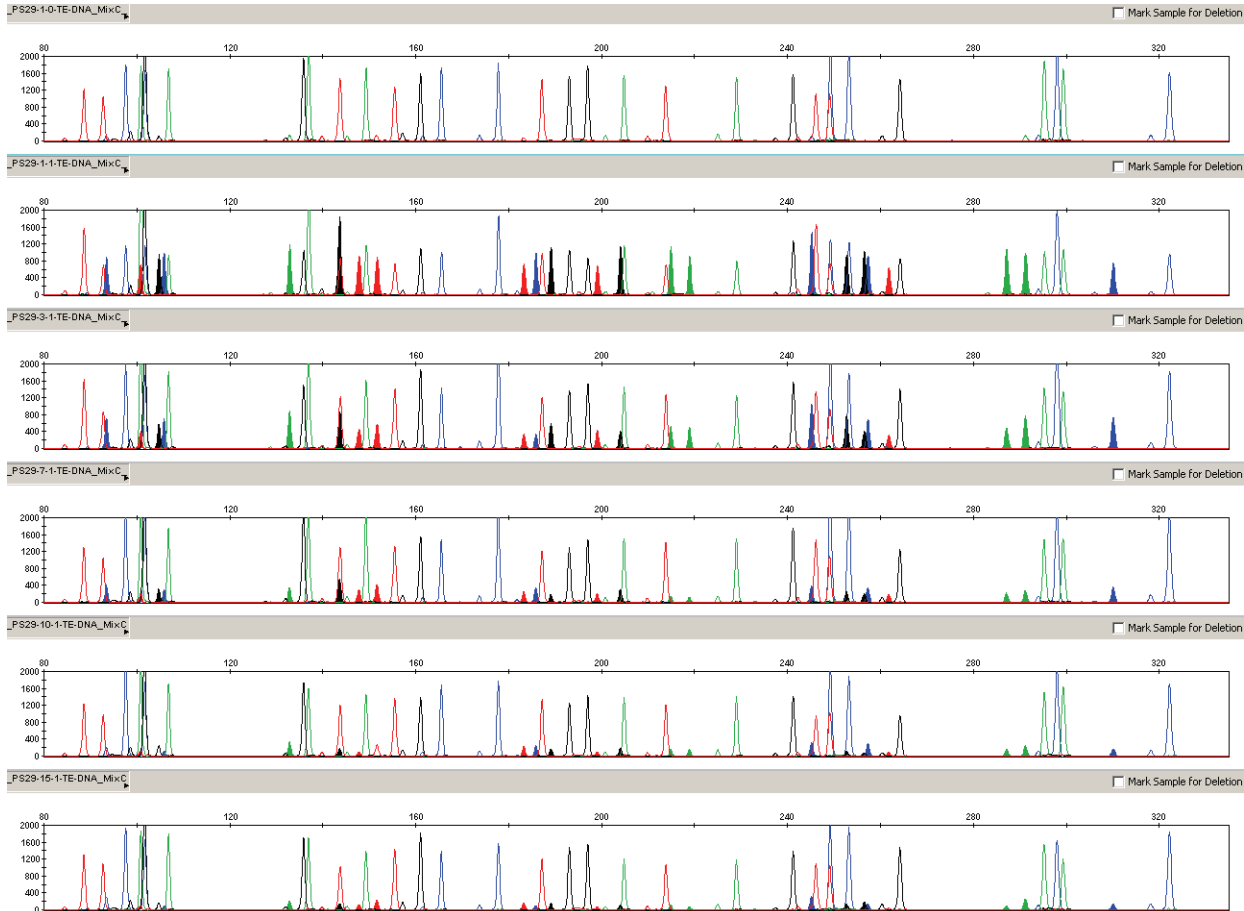


Table 9 Genotypes of mixed DNA samples

Locus	Sample A Genotype	Sample B Genotype
D10S1248	13, 14	12, 15
vWA	14, 17	17, 19
D16S539	10, 11	9, 12
D2S1338	17, 23	17, 20
AMEL	X, Y	X
DS1179	11,14	10,11
D21S11	29, 35	31.2, 32.2

Locus	Sample A Genotype	Sample B Genotype
D18S51	15, 16	13, 14
D22S10	45, 15	15, 16
D19S433	11, 17.2	13
TH01	7, 8	6, 9.3
FGA	19, 25	22, 23
D2S441	11, 12	11, 14
D3S1358	14, 17	15, 16
D1S1656	12, 18.3	11, 15
D12S391	18, 18.3	18, 22

Population data

SWGDM guideline 2.7

“The distribution of genetic markers in populations should be determined in relevant population groups.” (SWGDM, July 2003)

Overview

The NGM™ Kit contains loci for which extensive population data are available. For additional information on 11 loci shared between the kits, see the population data and additional studies section of the AmpF ℓ STR™ SGM Plus™ PCR Amplification Kit User's Manual (Part no. 44309589).

Population samples used in these studies

The NGM™ Kit was used to generate the population data provided in this section. Whole blood samples, provided by the Interstate Blood Bank (Memphis, Tennessee), were collected in the United States (with no geographical preference) from randomly selected individuals of known ethnicities. Ethnicities of sample donors were:

- African-American – 344 samples
- Caucasian – 346 samples
- Hispanic – 390 samples

DNA was extracted using an ABI PRISM™ 6100 Nucleic Acid PrepStation.

In addition to the alleles that were observed and recorded in our databases, other alleles have been published or reported to us by other laboratories (see the STRBase at www.cstl.nist.gov/div831/strbase).

New Primers added to the NGM™ Kit

Comparison of primer sequences with other AmpF ℓ STR™ Kits

Both the NGM™ and NGM SElect™ Kits contain an additional unlabeled primer for the D8S1179 locus to allow detection of a rare population-specific SNP-containing allele affecting one of the primer binding regions for that locus. The additional D8 primer was first introduced in the Identifiler™ kit (released 2001) and has been included in all other AmpF ℓ STR™ kits containing the D8S1179 locus since that time. This primer is not included in AmpF ℓ STR™ kits released before 2001, including the SGM Plus™ kit. Laboratories may therefore see occasional non-concordance at the D8S1179 locus when comparing results from different AmpF ℓ STR™ kits.

Both the NGM™ and NGM SElect™ Kits also contain slight modifications to the Amelogenin primers to reduce inter-species cross-reactivity compared to previous AmpFℓSTR™ kits.

Inclusion of three SNP-specific primers to address mutations at the Amelogenin, D2S441, and D22S1045 loci

After the initial release of the NGM™ Kit in early 2010, more recent population study results showed the existence of certain mutations that affected primer binding sites at three of the NGM™ Kit loci: amelogenin, D2S441, and D22S1045 (Carolyn Hill and John Butler, National Institute of Standards and Technology, personal communication). The mutations, when present, caused the drop-out of affected alleles. While the mutations are relatively rare and restricted primarily to specific population groups, the decision was made to include one additional PCR primer for each of the affected loci to allow the mutant alleles to be detected by the NGM™ Kit. The inclusion of the additional primers also required a minor re-optimization of the NGM™ Kit Master Mix to support the expanded primer mix.

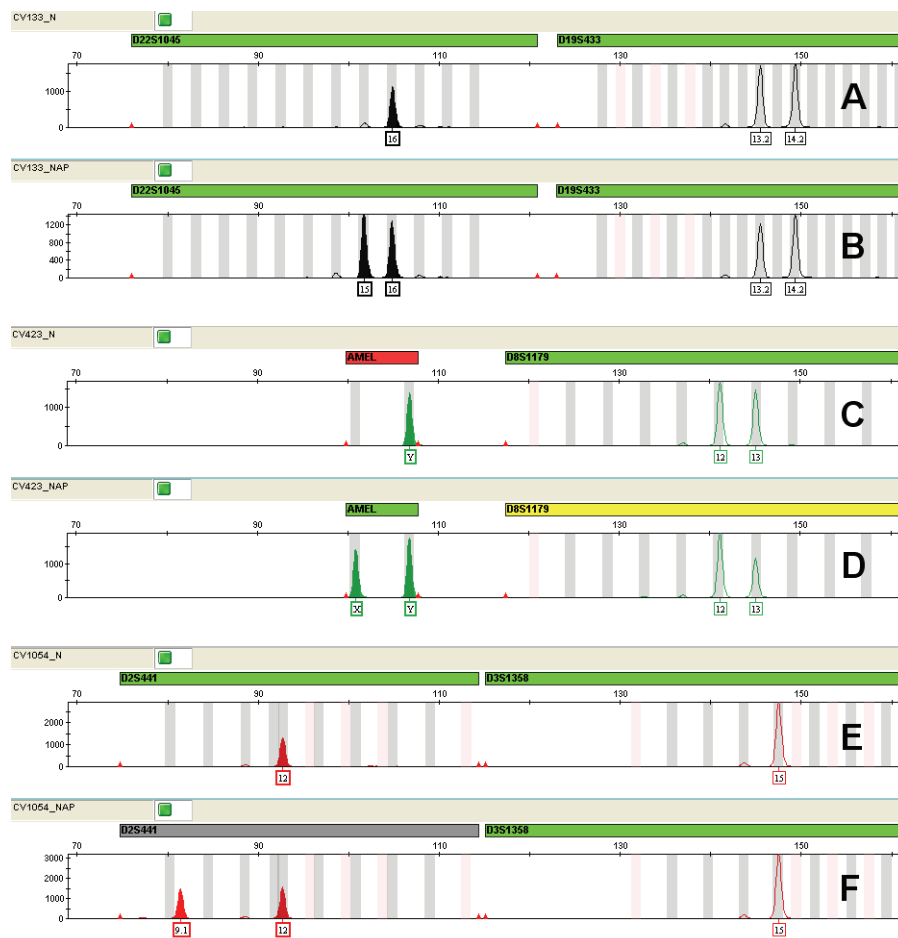
Subsequent validation experiments showed that other aspects of kit performance remained unaffected by the additional primers and the minor re-optimization of the Master Mix. Current versions of both the AmpFℓSTR™ NGM™ and NGM SElect™ Kits contain identical sets of primers for the 16 loci they have in common; the only difference between the kits is the presence of primers for the SE33 locus in the NGM SElect™ Kit.

The first commercially available batches of the NGM™ Kit to include the additional primers and benefit from the minor Master Mix re-optimisation were 200 Reaction Kit Lot 1105011 (Part no. 4415020) and 1000 Reaction Kit Lot 1106009 (Part no. 4415021). These lots were released in late summer 2011.

Note: The inclusion of the three additional primers in the NGM SElect™ Kit and the associated buffer optimisation studies were conducted as part of the original development of the NGM SElect™ Kit.

Figure 25 shows in the new version of the NGM™ Kit electropherograms with examples where the additional primers for amelogenin, D2S441 and, D22S1045 allow mutant alleles at these loci to be detected, whereas they had not been detected by the original version of the NGM™ Kit.

Figure 25 New PCR primers added to D22S1045, amelogenin and D2S441 loci. The electropherograms below show the effects of adding additional PCR primers to amplify known mutant alleles that would otherwise not be detected. Panels A and B show the D22S1045 locus for individual CV133 without (A) and with (B) the extra primer; the additional primer allowed the variant of allele 15 to be detected. Panels C and D show the amelogenin locus for individual CV423 without (C) and with (D) the extra primer; the additional primer allowed the variant of the X allele to be detected. Panels E and F show the D2S441 locus for individual CV1054 without (E) and with (F) the extra primer; when amplified, the variant allele was always seen to type as a 9.1 microvariant



Mutation rate

Estimating germ-line mutations

Estimation of spontaneous or induced germ-line 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.

In previous studies, genotypes of ten STR loci that were amplified by the AmpF Φ STRTM SGM PlusTM PCR Amplification 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; mutations were not detected at any of the other nine STR loci. The D18S11 mutation was represented by an increase of one 4-nt repeat unit, allele 17 was inherited as allele 18 (single-step mutation). The maternal/paternal source of this mutation could not be distinguished.

**Additional
mutation studies**

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 NGM™ 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 10 Allele frequencies (%) by population group for NGM™ Kit STR loci. (The + symbol indicates alleles not detected or, where values appear in parentheses, alleles not detected in significant quantities.)

Allele	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
D10S1248			
6	(0.15) [†]	+	+
7	(0.15) [†]	+	+
8	+	+	(0.13) [†]
9	(0.15) [†]	+	(0.13) [†]
10	(0.15) [†]	+	(0.13) [†]
11	3.63	(0.58) [†]	(0.26) [†]
12	13.95	3.47	4.23
13	22.67	29.05	25.51
14	28.2	29.77	36.03
15	18.6	19.65	22.95
16	9.88	13.44	8.08
17	2.18	3.76	2.56
18	(0.29) [†]	(0.29) [†]	+
19	+	+	+
20	+	+	+
D12S391			
13	+	+	(0.13) [†]
14	+	+	(0.13) [†]
15	7.12	4.19	3.97
15.1	(0.15) [†]	+	+
16	5.09	3.47	5.13
16.1	(0.15) [†]	+	+
17	15.7	10.55	7.31
17.1	(0.44) [†]	+	(0.26) [†]

Allele <i>(continued)</i>	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
17.3	(0.58) [†]	2.02	1.15
18	25.73	16.18	20
18.3	1.16	2.17	2.05
19	14.1	12.28	18.59
19.1	(0.58) [†]	†	†
19.3	(0.29) [†]	(0.58) [†]	1.15
20	11.92	9.83	17.31
20.3	†	†	†
21	6.98	13.73	8.72
21.3	(0.15) [†]	(0.14) [†]	†
22	5.09	10.69	6.92
23	3.34	8.09	3.72
24	(0.58) [†]	3.61	1.79
25	0.87	2.02	1.28
26	†	(0.29) [†]	(0.26) [†]
27	†	(0.14) [†]	(0.13) [†]
28	†	†	†
D16S539			
5	†	†	†
6	†	†	(0.13) [†]
7	†	†	†
8	3.49	1.45	1.92
9	22.24	12.43	10.38
10	11.63	4.05	15.77
11	29.07	32.23	31.92
12	19.33	30.78	24.49
12.2	†	†	†
13	12.94	17.2	13.97
14	1.31	1.73	1.15
15	†	(0.14) [†]	(0.26) [†]
16	†	†	†
D18S51			
7	†	†	†
8	†	†	†
9	†	†	(0.13) [†]
9.2	†	†	†
10	(0.29) [†]	1.16	0.64

Allele <i>(continued)</i>	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
10.2	(0.15) [†]	†	†
11	(0.44) [†]	0.87	1.15
11.2	†	†	†
12	6.25	14.74	10.77
12.2	†	†	†
13	4.07	11.85	11.54
13.2	(0.29) [†]	†	†
14	5.81	17.49	15.77
14.2	(0.44) [†]	†	(0.13) [†]
15	17.3	15.32	12.31
15.2	†	†	(0.13) [†]
16	18.31	11.85	12.82
16.2	†	†	†
17	15.7	10.98	17.05
17.2	†	†	†
18	14.1	8.53	8.46
18.2	†	†	†
19	9.16	4.34	3.46
19.2	†	†	†
20	4.22	1.3	1.79
20.2	(0.15) [†]	†	†
21	2.03	1.01	2.31
21.2	†	†	†
22	1.02	(0.29) [†]	0.64
22.2	†	†	†
23	(0.29) [†]	(0.29) [†]	(0.51) [†]
23.2	†	†	†
24	†	†	(0.26) [†]
25	†	†	(0.13) [†]
26	†	†	†
27	†	†	†
D19S433			
9	(0.29) [†]	†	†
9.2	†	†	†
10	1.16	(0.14) [†]	(0.38) [†]
10.2	(0.15) [†]	†	†
11	9.74	†	1.54

Allele <i>(continued)</i>	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
11.2	(0.29) [†]	†	(0.26) [†]
12	10.76	7.23	8.46
12.1	†	(0.14) [†]	†
12.2	3.78	(0.14) [†]	1.41
13	28.63	27.46	18.72
13.2	5.23	1.59	7.05
14	18.31	34.68	30.38
14.2	5.81	2.17	4.49
15	6.25	16.18	12.69
15.2	4.22	3.47	7.31
16	1.45	5.92	4.1
16.2	3.34	(0.29) [†]	2.31
17	†	(0.29) [†]	(0.51) [†]
17.2	(0.58) [†]	(0.14) [†]	(0.38) [†]
18	†	(0.14) [†]	†
18.2	†	†	†
D1S1656			
9	(0.15) [†]	†	(0.13) [†]
10	1.31	(0.29) [†]	(0.38) [†]
11	5.52	5.92	3.85
12	7.99	16.04	9.49
13	11.19	6.65	7.18
14	24.27	6.36	11.03
14.3	1.02	(0.29) [†]	(0.26) [†]
15	17.44	15.17	16.03
15.3	1.89	8.53	2.95
16	10.32	9.97	15.26
16.1	†	†	(0.26) [†]
16.3	7.27	4.91	5.26
17	2.62	4.91	6.92
17.1	†	(0.29) [†]	†
17.3	5.96	12.72	15.26
18	(0.44) [†]	(0.29) [†]	0.77
18.3	1.89	5.92	4.36
19	(0.15) [†]	†	†
19.3	(0.58) [†]	1.73	0.64
20.3	†	†	†

Allele (continued)	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
D21S11			
23.2	†	†	†
24	†	†	†
24.2	†	†	-- (0.26)
25	†	†	†
25.2	†	†	†
26	(0.29) [†]	(0.58) [†]	(0.38) [†]
26.2	†	†	†
27	5.52	2.75	1.41
27.1	(0.15) [†]	†	†
27.2	†	†	†
28	25.29	16.91	11.03
28.2	†	†	(0.13) [†]
29	15.7	23.55	21.15
29.2	†	(0.14) [†]	†
29.3	(0.15) [†]	(0.14) [†]	†
30	20.93	23.41	27.95
30.2	1.6	2.75	1.67
31	8.72	6.79	5
31.2	4.94	8.67	11.28
32	1.31	2.31	1.28
32.2	6.98	9.54	12.44
33	0.87	†	-- (0.13)
33.2	3.2	2.02	5.26
34	(0.15) [†]	†	†
34.2	†	(0.43) [†]	(0.13) [†]
35	3.49	†	(0.26) [†]
35.2	†	†	†
36	0.73	†	(0.13) [†]
36.2	†	†	†
37	†	†	†
37.2	†	†	†
38	†	†	(0.13) [†]
38.2	†	†	†
39	†	†	†
D22S1045			
5	†	†	†

Allele (continued)	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
6	†	†	†
7	†	†	†
8	0.73	†	†
9	†	†	†
10	4.07	(0.43) [†]	0.64
11	14.53	13.58	7.82
12	5.96	(0.58) [†]	1.03
13	(0.29) [†]	1.01	1.03
14	8.58	3.47	2.18
15	23.55	36.56	42.56
16	19.91	36.27	35.64
17	20.35	7.51	7.95
18	2.03	(0.58) [†]	1.03
20	†	†	(0.13) [†]
D2S1338			
13	(0.15) [†]	(0.14) [†]	†
14	†	(0.14) [†]	†
15	(0.29) [†]	(0.14) [†]	†
16	5.23	4.19	3.59
17	10.03	18.79	17.69
18	4.8	8.38	6.54
19	15.99	14.31	17.82
20	10.03	15.46	13.72
21	12.79	2.75	3.59
22	12.65	1.73	6.28
23	9.3	10.12	14.87
24	8.58	9.97	8.85
25	6.98	11.85	5.38
26	2.47	1.73	1.41
27	0.73	(0.29) [†]	(0.13) [†]
28	†	†	(0.13) [†]
29	†	†	†
D2S441			
8	(0.15) [†]	†	†
9	†	(0.58) [†]	(0.26) [†]
10	8.87	19.8	31.15
11	35.03	33.82	31.67

Allele <i>(continued)</i>	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
11.3	3.34	5.06	4.36
12	20.06	4.05	3.97
12.3	(0.15) [†]	(0.29) [†]	(0.38) [†]
13	3.63	3.18	1.92
13.3	†	†	†
14	26.89	28.32	22.56
14.3	†	†	†
15	1.89	4.48	3.33
16	†	(0.26) [†]	(0.38) [†]
17	†	†	†
D3S1358			
9	(0.29) [†]	†	(0.13) [†]
11	†	(0.29) [†]	†
12	(0.15) [†]	†	(0.13) [†]
13	(0.58) [†]	(0.14) [†]	(0.38) [†]
14	9.16	15.17	9.49
15	28.34	27.31	34.49
15.2	(0.29) [†]	†	†
16	32.85	23.99	26.54
16.2	†	†	†
17	22.09	19.8	17.95
17.2	†	†	†
18	5.81	11.85	10.13
18.2	†	†	†
19	(0.44) [†]	1.45	0.77
20	†	†	†
D8S1179			
7	†	†	†
8	(0.44) [†]	2.02	0.64
9	(0.29) [†]	1.3	(0.26) [†]
10	3.34	10.84	9.49
11	5.81	6.65	4.87
12	11.05	15.03	12.44
13	18.31	33.53	33.33
14	36.05	18.64	23.46
15	17.73	8.67	11.54
16	5.96	2.89	3.33

Allele <i>(continued)</i>	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
17	1.02	(0.43) [†]	0.64
18	†	†	†
19	†	†	†
20	†	†	†
FGA			
16	†	(0.14) [†]	†
16.1	(0.29) [†]	†	†
16.2	†	†	†
17	†	(0.14) [†]	†
17.2	†	†	†
18	0.87	1.16	0.64
18.2	0.73	†	†
19	6.83	5.35	7.82
19.2	(0.44) [†]	†	†
20	6.69	15.61	8.72
20.2	(0.29) [†]	(0.43) [†]	(0.26) [†]
21	12.06	18.35	13.59
21.2	(0.15) [†]	(0.29) [†]	†
22	18.17	18.93	14.1
22.2	(0.15) [†]	0.87	0.64
23	16.86	14.6	12.69
23.2	†	(0.43) [†]	(0.38) [†]
23.3	(0.29) [†]	†	†
24	18.75	14.31	15.9
24.2	†	†	†
25	9.3	6.79	14.1
25.2	†	†	†
26	3.92	1.88	6.79
26.2	†	†	†
27	2.62	(0.58) [†]	2.95
27.2	†	†	†
28	1.16	(0.14) [†]	0.9
--	†	†	†
29	†	†	(0.38) [†]
29.2	†	†	†
30	(0.15) [†]	†	(0.13) [†]
30.2	(0.15) [†]	†	†

Allele <i>(continued)</i>	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
31	†	†	†
31.2	†	†	†
32	†	†	†
32.2	†	†	†
33.2	†	†	†
34.2	(0.15) [†]	†	†
42.2	†	†	†
43.2	†	†	†
44.2	†	†	†
45.2	†	†	†
46.2	†	†	†
47.2	†	†	†
48.2	†	†	†
49.2	†	†	†
50.2	†	†	†
51.2	†	†	†
TH01			
3	†	†	†
4	†	†	†
5	(0.44) [†]	(0.14) [†]	†
5.3	†	†	†
6	15.41	21.68	27.95
6.1	(0.15) [†]	†	†
6.3	†	†	†
7	37.06	17.77	31.92
7.3	†	†	†
8	21.22	11.42	8.46
8.3	†	†	†
9	15.84	17.05	12.56
9.3	8.43	31.07	17.69
10	1.45	0.87	1.41
10.3	†	†	†
11	†	†	†
12	†	†	†
13	†	†	†
13.3	†	†	†

Allele (continued)	African American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
vWA			
10	†	†	†
11	(0.44) [†]	†	(0.13) [†]
12	†	†	(0.26) [†]
13	0.73	(0.14) [†]	(0.13) [†]
14	7.41	8.38	6.41
15	20.93	12.14	9.87
15.2	†	†	†
16	27.47	22.69	30
17	19.48	27.31	27.18
17.3	†	†	(0.13) [†]
18	13.81	18.06	18.46
18.2	†	†	†
19	6.98	9.97	6.67
20	2.03	1.3	0.77
21	(0.58) [†]	†	†
22	†	†	†
23	(0.15) [†]	†	†
24	†	†	†
25	†	†	†

† A minimum allele frequency (0.7% for the African-American database, 0.7% for the U.S. Caucasian database, 0.9% for the U.S. Hispanic database, and 1.3% for the Native American database) is suggested by the National Research Council in forensic calculations.

Table 11 shows the allele frequencies at NGM™ Kit loci by population group. The P_I value is the probability that two individuals selected at random will have an identical NGM™ Kit genotype (Sensabaugh, 1982). The P_I values for the populations described in this section are then 6.74 × 10⁻²⁰ ng/μL⁻²² (African-American), 2.76 × 10⁻¹⁹ (U.S. Caucasian) and 4.00 × 10⁻¹⁹ (U.S. Hispanic).

Table 11 Probability of identity (P_I) values for the NGM™ Kit STR loci

Locus	African-American (N = 344)	U.S.-Caucasian (N = 346)	U.S.-Hispanic (N = 390)
D10S1248	0.069	0.094	0.111
vWA	0.062	0.065	0.091
D16S539	0.072	0.104	0.082
D2S1338	0.023	0.032	0.032
D8S1179	0.076	0.063	0.068
D21S11	0.045	0.052	0.050

Locus	African-American (N = 344)	U.S.-Caucasian (N = 346)	U.S.-Hispanic (N = 390)
D18S51	0.031	0.031	0.028
D22S1045	0.056	0.133	0.161
D19S433	0.040	0.085	0.048
TH01	0.094	0.080	0.091
FGA	0.033	0.039	0.028
D2S441	0.101	0.098	0.107
D3S1358	0.100	0.075	0.095
D1S1656	0.034	0.022	0.025
D12S391	0.039	0.023	0.032
Combined	6.74×10^{-20}	2.76×10^{-19}	4.00×10^{-19}

Probability of paternity exclusion

Table 12 Probability of Paternity Exclusion values for the NGM™ Kit™ STR loci

Locus	African-American (N = 344)	Caucasian (N = 346)	Hispanic (N = 390)
D10S1248	0.659	0.568	0.469
vWA	0.603	0.644	0.638
D16S539	0.556	0.579	0.557
D2S1338	0.798	0.752	0.738
D8S1179	0.587	0.616	0.557
D21S11	0.727	0.729	0.692
D18S51	0.786	0.758	0.718
D22S1045	0.698	0.455	0.397
D19S433	0.636	0.507	0.647
TH01	0.515	0.502	0.539
FGA	0.727	0.666	0.708
D2S441	0.522	0.497	0.477
D3S1358	0.476	0.522	0.482
D1S1656	0.745	0.811	0.723
D12S391	0.745	0.799	0.662
Combined	0.999999931	0.999999835	0.999999376

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 NGM™ Kit STR loci (Chakraborty, Stivers, and Zhong, 1996).



Troubleshooting

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

Observation	Possible causes	Recommended actions
Faint or no signal from both, the 007 AmpF ζ STR TM Control DNA and the DNA test samples at all loci	Incorrect volume or absence of either AmpF ζ STR TM NGM TM Master Mix or AmpF ζ STR TM NGM TM Primer Set	Repeat amplification using correct reagent volumes.
	No activation of enzyme	Repeat amplification, making sure to hold reactions initially at 95°C for 11 min.
	Master Mix not vortexed thoroughly before aliquoting	Vortex Master Mix thoroughly.
	AmpF ζ STR TM NGM TM Primer Set exposed to too much light	Store Primer Set protected from light.
	GeneAmp TM PCR System malfunction	Refer to the thermal cycler user's manual and check instrument calibration.
	Incorrect thermal cycler parameters	Check the protocol for correct thermal cycler parameters.
	Tubes/plate not seated tightly in the thermal cycler during amplification	Push reaction tubes/plate firmly into contact with block after first cycle. Repeat test.
	Wrong PCR reaction tubes or plate	Use Applied Biosystems TM MicroAmp Reaction Tubes with Caps or the MicroAmp Optical 96-Well Reaction Plate for the GeneAmp TM PCR System 9700 or Veriti TM 96-well Thermal Cycler.
	MicroAmp TM Base used with tray/retainer set and tubes in GeneAmp TM PCR System 9700	Remove MicroAmp Base from tray/retainer set and repeat test.
	Insufficient PCR product electrokinetically injected	Refer to Chapter 3, "Electrophoresis" on page 25, for instructions on recommended actions on the ABI PRISM TM 3100/ 3100- <i>Avant</i> or Applied Biosystems TM 3130/3130 <i>xl</i> , 3500/ 3500 <i>xl</i> , and the ABI PRISM TM 310 instruments.
Degraded formamide	Check the storage of formamide; do not thaw and refreeze multiple times. Try Hi-Di TM Formamide.	

Observation	Possible causes	Recommended actions
Positive signal from AmpF \mathcal{L} STR™ Control DNA 007 but partial or no signal from DNA test samples	Quantity of test DNA sample is below assay sensitivity	Quantify DNA and add either 1.0 ng or 500 pg of DNA depending upon cycle number being used. Repeat test.
	Test sample contains high concentration of PCR inhibitor (for example, heme compounds, certain dyes)	Quantify 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 \mathcal{L} 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 two alleles present at a locus	Presence of exogenous DNA	Use appropriate techniques to avoid introducing foreign DNA during laboratory handling.
	Amplification of stutter product	Interpret according to laboratory procedures. Note: Additional information will be provided on completion of validation.
	Mixed sample	
	Incomplete 3' A base addition (n-1 nt position)	Addition of excess DNA to the reaction will contribute to the occurrence of incomplete 3' base addition. Quantify DNA and add 1.0 ng of DNA to the reaction. Repeat test. Also be sure to include the final extension step of 60°C for 10 min in the PCR.
	Signal exceeds dynamic range of instrument (off-scale data)	Ensure cycle number is optimized according to instructions on page 22. Repeat PCR amplification using fewer PCR cycles or use your laboratory's SOP to analyze off-scale data.
	Poor spectral separation (bad matrix)	Follow the steps for creating a spectral file.
		Confirm that Filter Set G5 modules are installed and used for analysis.
	Too much DNA in reaction	Use recommended amount of template DNA: 1.0 ng at 29 cycles; 500 pg at 30 cycles.
Incomplete denaturation of double stranded DNA	Use the recommended amount of Hi-Di™ Formamide and perform heat denaturation according to instructions in Chapter 3, "Electrophoresis".	
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, gold-plated silver blocks or Veriti™ 96-well Thermal Cycler only.



Ordering Information

Materials and equipment not included

The tables below list optional equipment and materials not supplied with the NGM™ Kit. Unless otherwise noted, many of the items are available from major laboratory suppliers (MLS).

Equipment	Part number
Applied Biosystems™ 3500/3500xL Genetic Analyzer for Human Identification	Contact your local Life Technologies sales representative
ABI PRISM™ 3100/3100- <i>Avant</i> Genetic Analyzer	
Applied Biosystems™ 3130/3130 <i>xl</i> Genetic Analyzer	
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
Silver 96-Well Sample Block	N8050251
Gold-plated Silver 96-Well Sample Block	4314443
Veriti™ 96-well Thermal Cycler	4375786
ProFlex™ 96-Well PCR System	4484075
Tabletop centrifuge with 96-Well Plate Adapters (optional)	MLS

Item	Part number
3500/3500xL Analyzer materials	
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™ 600 LIZ™ Size Standard v2.0	4408399

Item	Part number
Note: 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> (Part no. 4401661)	
AmpF _{STR} [™] NGM [™] PCR Amplification Kit (200x/1000x)	4415020/4415021
3100/3100-Avant Analyzer materials	
96-Well Plate Septa	4315933
Reservoir Septa	4315932
3100/3100-Avant Genetic Analyzer Capillary Array, 36-cm	4333464
POP-4 [™] Polymer for 3100/3100-Avant Genetic Analyzers	4316355
3100/3100-Avant Genetic Analyzer Autosampler Plate Kit, 96-well	4316471
GeneScan [™] 500 LIZ [™] Size Standard <i>Or</i>	4322682
GeneScan [™] 600 LIZ [™] Size Standard v2.0	<i>Or</i> 4408399
Running Buffer, 10X	402824
DS-33 Matrix Standard Kit (Dye Set G5)	4345833
MicroAmp [™] Optical 96-Well Reaction Plate	N8010560
250-µL Glass Syringe (array-fill syringe)	4304470
5.0-mL Glass Syringe (polymer-reserve syringe)	628-3731
Note: For a complete list of parts and accessories for the 3100 instrument, refer to Appendix B of the <i>ABI PRISM™ 3100 Genetic Analyzer and 3100-Avant Genetic Analyzer User Reference Guide</i> (Part no. 4335393).	
3130/3130xl Analyzer materials	
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
3130/3130xl Genetic Analyzer Autosampler Plate Kit, 96-well	4316471
GeneScan [™] 500 LIZ [™] Size Standard <i>Or</i>	4322682
GeneScan [™] 600 LIZ [™] Size Standard v2.0	<i>Or</i> 4408399
Running Buffer, 10X	402824
DS-33 Matrix Standard Kit (Dye Set G5)	4345833
MicroAmp [™] Optical 96-Well Reaction Plate	N8010560
For a complete list of parts and accessories for the 3130xl instrument, refer to Appendix A of the <i>Applied Biosystems 3130/3130xl Genetic Analyzers Maintenance, Troubleshooting, and Reference Guide</i> (Part no. 4352716).	
310 DNA Analyzer materials	
310 DNA Analyzer Capillary Array, 47-cm	402839
0.5 mL Sample Tray	5572
96-Well Tray Adaptor (for 9700 thermal cycler trays)	4305051

Item	Part number
GeneScan™ 500 LIZ™ Size Standard <i>Or</i>	4322682
GeneScan™ 600 LIZ™ Size Standard v2.0	<i>Or</i> 4408399
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-33 Matrix Standard Set (6-FAM™, VIC™, NED™, PET™, and LIZ™ dyes) for ABI PRISM™ 310/377 systems	4318159
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>ABI PRISM™ 310 Genetic Analyzer User Guide</i> (Part 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

B

Appendix B Ordering Information
Materials and equipment not included



PCR Work Areas

- Work area setup and lab design 111
- PCR setup work area 111
- Amplified DNA work area 112

Work area setup and lab design

Many resources are available for the appropriate design of a PCR laboratory. If you are using the AmpF ϕ STR™ NGM™ PCR Amplification Kit for:

- Forensic DNA testing, refer to “Forensic Laboratories: Handbook for Facility Planning, Design, Construction and Moving,” National Institute of Justice, 1998 (www.nij.org/publications)
- 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 NGM™ Kit (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 NGM™ 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 NGM™ Kit.

- Veriti™ 96-well Thermal Cycler



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.
-

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
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030), found at: 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

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/





Documentation and Support

Related documentation

The following related documents are shipped with the system:

Document	Part number
<i>ABI PRISM™ 3100/3100-Avant Data Collection v2.0 User Guide</i>	4347102
<i>ABI PRISM™ 3100/3100-Avant Genetic Analyzers Using Data Collection Software v2.0 User Bulletin</i>	4350218
<i>ABI PRISM™ 3100 Genetic Analyzer User Manual (Data Collection v1.1)</i>	4315834
<i>ABI PRISM™ 3100/3100-Avant Genetic Analyzers Protocols for Processing AmpFtSTR™ PCR Amplification Kit PCR Products User Bulletin</i>	4332345
<i>AmpFtSTR™ NGM™ PCR Amplification Kit - PCR Setup Quick Reference Card</i>	4442401
<i>AmpFtSTR™ NGM™ PCR Amplification Kit - CE Quick Reference Card</i>	4442693
<i>Veriti™ 96-Well Thermal Cycler AmpFtSTR™ Kit Validation User Bulletin</i>	4440754
<i>ProFlex™ PCR System Kit Validation User Bulletin</i>	100031595
<i>Applied Biosystems™ 3130/3130xl 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™ 3130/3130xl DNA Analyzers User Guide</i>	4331468
<i>Applied Biosystems™ 3730/3730xl Genetic Analyzer Getting Started Guide</i>	4359476
<i>Quantifiler™ Kits: Quantifiler™ Human DNA Quantification Kit and Quantifiler™ Y Human Male DNA Quantification Kit User's Manual</i>	4344790
<i>Identifiler™ 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

Document	Part number
<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
<i>GeneMapper™ ID-X Software Version 1.1 (Mixture Analysis) Quick Reference Guide</i>	4402094
<i>GeneMapper™ ID-X Software Version 1.2 Quick Reference Guide</i>	4426482
<i>GeneMapper™ ID-X Software Version 1.2 Reference Guide</i>	4426481

Note: To open the user documentation, use the Adobe™ Reader™ software available from www.adobe.com

Note: For additional documentation, see “Obtain support” on page 118.

Obtain support

For HID support:

- **In North America** – Send an email to 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

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 their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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.
- American Association of Blood Banks. 2004. *Guidance for Standards for Parentage Relationship Testing Laboratories*. 7th ed. Bethesda, Md: American Association of Blood Banks. 58 pp.
- 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. *Int. J. Leg. Med.* 108: 31-35.
- 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.
- 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. *Int. J. Leg. Med.* 108: 180-185.
- Baron, H., Fung, S., Aydin, A., Bahrig, S., Luft, F.C., Schuster, H. 1996. Oligonucleotide ligation assay (OLA) for the diagnosis of familial hypercholesterolemia. *Nat. Biotechnol.* 14:1279–1282.
- Begovich A.B., McClure G.R., Suraj V.C., Helmuth R.C., Fildes N., Bugawan T.L., Erlich H.A., Klitz W. 1992. Polymorphism, recombination, and linkage disequilibrium within the HLA class II region. *J. Immunol.* 148:249–58.
- Bender, K., Farfan, M.J., Schneider, P.M. 2004. Preparation of degraded human DNA under controlled conditions. *Forensic Sci. Int.* 139:134–140.
- Bonferroni, C.E. 1936. Teoria statistica delle classi e calcolo delle probabilita. *Publicazioni del R Istituto Superiore di Scienze Economiche e Commerciali di Firenze* 8:3–62.
- Brinkman, B., Klintschar, 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.
- Brinkman, B., Moller, A. and Wiegand, P. 1995. Structure of new mutations in 2 STR systems. *Intl. J. Legal Med.* 107:201–203.
- Butler, J.M. 2005. *Forensic DNA Typing*. Burlington, MA:Elsevier Academic Press.
- Butler, J.M., Shen, Y., McCord, B.R. 2003. The development of reduced size STR amplicons as tools for analysis of degraded DNA. *J. Forensic Sci.* 48:1054–1064.
- 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.

- 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. and Stivers, D.N. 1996. Paternity exclusion by DNA markers: effects of paternal mutations. *J. Forensic Sci.* 41:671–677.
- Chung, D.T., Drabek, J., Opel, K.L., Butler, J.M. and McCord, B.R. 2004. A study of the effects of degradation and template concentration on the amplification efficiency of the Miniplex primer sets. *J. Forensic Sci.* 49:733–740.
- Clark J.M. 1988. Novel non-templated nucleotide addition reactions catalyzed by procaryotic and eucaryotic DNA polymerases. *Nucleic Acids Res.* 16:9677–9686.
- Coble, M.D. and Butler, J.M. 2005. Characterization of new miniSTR loci to aid analysis of degraded DNA. *J. Forensic Sci.* 50:43–53.
- 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.
- Drabek, J., Chung, D.T., Butler, J.M., McCord, B.R. 2004. Concordance study between Miniplex assays and a commercial STR typing kit. *J. Forensic Sci.* 49:859–860.
- 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.
- Farfan, M. J., Sanz, P., Lareu, M. V., and Carracedo, A. 9-30-1999. Population data on the D1S1656 and D12S391 STR loci in Andalusia (south Spain) and the maghreb (north Africa). *Forensic Sci.Int.* 104(1): 33-36.
- Federal Bureau of Investigation DNA Advisory Board. 1998. *Quality Assurance Standards for Forensic DNA Testing Laboratories*. Washington, DC: Federal Bureau of Investigation.
- Frank, W., Llewellyn, B., Fish, P., et al. 2001. Validation of the AmpF Φ STRTM Profiler PlusTM PCR Amplification Kit for use in forensic casework. *J. Forensic Sci.* 46:642–646.
- Gamero, J. J., Romero, J. L., Gonzalez, J. L., Arufe, M. I., Cuesta, M. I., Corte-Real, F., Carvalho, M., Anjos, M. J., Vieira, D. N., and Vide, M. C. 6-5-2000. A study on ten short tandem repeat systems: African immigrant and Spanish population data. *Forensic Sci.Int.* 110(3): 167-177.
- Gill, P., Fereday, L., Morling, N., Schneider, P.M. (2006) New multiplexes for Europe—Amendments and clarification of strategic development. *Forensic Sci. Int.* 163 (2006) 155–157.
- Glock, B., Dauber, E.M., Schwartz, D.W., Mayr W.R. 1997. Additional variability at the D12S391 STR locus in an Austrian population sample: sequencing data and allele distribution. *Forensic Sci. Int.* 90:197–203.
- Grossman, P.D., Bloch, W., Brinson, E., Chang, C.C., Eggerding, F.A., Fung, S., Iovannisci, D.M., Woo, S., Winn-Deen, E.S. 1994. High-density multiplex detection of nucleic acid sequences: oligonucleotide ligation assay and sequence-coded separation. *Nucleic Acids Res.* 22:4527–4534.

- Grubwieser, P., Muhlmann, R., Berger, B., Niederstatter, H., Palvic, M., Parson, W. 2006. A new "mini-STR-multiplex" displaying reduced amplicon lengths for the analysis of degraded DNA. *Int. J. Legal Med.* 120:115–120.
- Guo S.W., and Thompson, E.A. 1992. Performing the exact test of Hardy-Weinberg proportion for multiple alleles. *Biometrics* 48:361–372.
- Guthmiller, J.M., Vargas, K.G., Srikantha, R., Schomberg, L.L., Weistroffer, P.L., McCray, P.B. and Tack, B.F. 2001. Susceptibilities of oral bacteria and yeast to mammalian cathelicidins. *Antimicrob. Agents Chemother.* 45:3216–3219.
- 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.
- Holt, C., Stauffer, C., Wallin, J., et al. 2000. Practical applications of genotypic Surveys for forensic STR testing. *Forensic Sci. Int.* 112:91–109.
- Kalinowski, S.T. 2006. HW-QuickCheck: an easy-to-use computer program for checking genotypes for agreement with Hardy-Weinberg expectations. *Molecular Ecology Notes* 6:974–979.
- Kimpton, C., Walton, A., and Gill, P. 1992. A further tetranucleotide repeat polymorphism in the vWF gene. *Hum. Mol. Genet.* 1:287.
- Kong, X., Murphy, K., Raj, T., He, C., White, P.S., Matisse, T.C. 2004. A combined linkage-physical map of the human genome. *Am. J. Hum. Genet.* 75:1143–1148.
- Lareu, M.V., Barral, S., Salas, A., Pestoni, C., and Carracedo, A. 1998. Sequence variation of a hypervariable short tandem repeat at the D1S1656 locus. *Int. J. Legal Med.* 111(5):244–247.
- Lareu, M.V., Pestoni, M.C., Barros, F., Salas, A., Carracedo, A. 1996. Sequence variation of a hypervariable short tandem repeat at the D12S391 locus. *Gene* 182:151–153.
- 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.
- 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.
- Magnuson, V.L., Ally, D.S., Nylund, S.J., Karanjawala, Z.E., Rayman, J.B., Knapp, J.I., Lowe, A.L., Ghosh, S., 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.
- Mansfield, E.S., Robertson, J.M., Vainer, M., Isenberg, A.R., Frazier, R.R., Ferguson, K., Chow, S., Harris, D.W., Barker, D.L., Gill, P.D., Budowle, B., McCord, B.R. 1998. Analysis of multiplexed short tandem repeat (STR) systems using capillary array electrophoresis. *Electrophoresis* 19:101–107.
- 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. and Brinkmann, B. 1994. Locus ACTBP2 (SE33): Sequencing data reveal considerable polymorphism. *Int. J. Leg. Med.* 106: 262-267

- Moller, A. and Brinkmann, B. 1995. PCR-VNTRs (PCR-Variable Number of Tandem Repeats) in forensic science. *Cellular & Molec. Bio.* 41(5):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.
- Moretti, T., Baumstark, A., Defenbaugh, D., Keys, K., Smerick, J., and Budowle, B. 2001. Validation of short tandem repeats (STRs) for forensic usage: Performance testing of fluorescent multiplex STR systems and analysis of authentic and simulated forensic samples. *J. Forensic Sci.* 46(3):647-660.
- Mulero, J.J., Chang, C.W., and Hennessy, L.K. 2006. Characterization of N+3 stutter product in the trinucleotide repeat locus DYS392. *J. Forensic Sci.* 51:826-830.
- Nakahori, Y., Takenaka, O., and Nakagome, Y. 1991. A human X-Y homologous region encodes amelogenin. *Genomics* 9:264-269.
- National Institute of Justice Office of Law Enforcement Standards. 1998. *Forensic Laboratories: Handbook for Facility Planning, Design, Construction and Moving*. Washington, DC: National Institute of Justice. 76 pp.
- Puers C, Hammond HA, Jin L, Caskey CT, Schumm JW., Identification of repeat sequence heterogeneity at the polymorphic short tandem repeat locus HUMTH01[AATG]_n and reassignment of alleles in population analysis by using a locus-specific allelic ladder. 1: *Am J. Hum. Genet.* 1993 Oct;53(4):953-8
- Raymond M. & Rousset F., 1995. GENEPOP (version 1.2): population genetics software for exact tests and ecumenicism. *J. Heredity* 86:248-249.
- Revised Validation Guidelines-Scientific Working Group on DNA Analysis Methods (SWGDM). *Forensic Sci. Communications* (July 2004) Volume 6 (3). Available at www.fbi.gov/hq/lab/fsc/current/standards/2004_03_standards02.htm
- Rousset, F. 2008. Genepop'007: A complete reimplement of the Genepop software for Windows and Linux. *Molecular Ecology Resources* 8:103-106.
- Sensabaugh, G.F. 1982. Biochemical markers of individuality. In: Saferstein, R., ed. *Forensic Science Handbook*. Prentice-Hall, Inc., New York, pp. 338-415.
- Sharma, V., and Litt, M. 1992. Tetranucleotide repeat polymorphism at the D21S11 locus. *Hum Mol. Genet.* 1:67.
- Shin, C.H., Jang, P., Hong, K.M., Paik, M.K. 2004. Allele frequencies of 10 STR loci in Koreans. *Forensic Sci. Int.* 140:133-135.
- 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.
- 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.
- Suido, H., Nakamura, M., Mashimo, P.A., Zambon, J.J., and Genco, R.J. 1986. Arylaminopeptidase activities of the oral bacteria. *J. Dent. Res.* 65:1335–1340.
- Szibor, R., Lautsch, S., Plate, I., Bender, K., Krause, D. 1998. Population genetic data of the STR HumD3S1358 in two regions of Germany. *Int J Legal Med.* 111(3):160-1.
- Waiyawuth, W., Zhang, L., Rittner, C., Schneider, P.M. 1998. Genetic analysis of the short tandem repeat system D12S391 in the German and three Asian populations. *Forensic Sci. Int.* 94:25–31.
- Wallin, J.M., Buoncristiani, M.R., Lazaruk, K.D., Fildes, N., Holt, C.L., Walsh, P.S. 1998. SWGDAM validation of the AmpF \mathcal{L} STR™ blue PCR amplification kit for forensic casework analysis. *J. Forensic Sci.* 43:854–870.
- Wallin, J.M., Holt, C.L., Lazaruk, K.D., Nguyen, T.H., Walsh, P.S. 2002. Constructing universal multiplex PCR systems for comparative genotyping. *J. Forensic Sci.* 47:52–65.
- Walsh, P.S., Fildes, N.J., 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). 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.
- Wiegand, P., Schneider, H.R., Schurenkamp, M., Kleiber, M., and Brinkmann, B. 1998. Tetranucleotide STR system D8S1132: sequencing data and population genetic comparisons. *Int. J. Legal Med.* 111(4):180-182
- Wiegand, P. and Kleiber, M. 2001. Less is more—length reduction of STR amplicons using redesigned primers. *Int. J. Legal Med.* 114:285–287.
- Budowle B, Ge J, Chakraborty R, Eisenberg AJ, Green R, Mulero J, Lagace R, Hennessy L. 2011. Population genetic analyses of the NGM STR loci. *Int J Legal Med.* 125:101-9. Epub 2010 Sep.

Index

Symbols

- .fsa sample files 34, 46
- +A nucleotide addition
 - defined 76
 - efficiency of 76

Numerics

- 3100/3100-Avant instruments 27

A

- accuracy and reproducibility 63
- alleles
 - off-ladder 64
- alleles,off-ladder 64
- allelic ladder
 - about 18
 - figure 13
 - precision results table 65
 - requirements for accurate genotyping 25
 - volume per reaction 28, 30, 32
- amplification
 - loci 12
 - using bloodstained FTA cards 23
- annealing temperatures, validation of 61, 62
- artifacts 77

B

- baseline noise, examples 78
- biohazard safety 114

C

- CEPH 80
- characterization of loci, validation 79
- chemical safety 114
- contents of kit 17, 20
- control DNA
 - about 18
- control DNA 007 14, 18
- cycle number, validation 62

D

- data
 - accuracy, precision, reproducibility 63
 - for different populations 91
- Data Collection Software 16
- developmental validation 60
- DNA
 - control, about 18
 - degraded 84
 - effect of quantity 83
 - mixture studies 88
 - mixture studies figure 89
 - mixtures, limit of detection 89
 - negative-control reaction 22
 - positive-control reaction 22
 - quantification methods 19, 20
 - sample preparation 22
 - test sample 22
- documentation, related 117

E

- electrophoresis
 - Data Collection Software 27, 29, 31
 - prepare samples 28
 - preparing samples on the 310 instrument 31
 - preparing samples on the 3500/3500xL instrument 29
 - reagents and parts 29, 31
 - references 27, 29, 31
 - run module 27, 29, 31
 - set up 29, 31
 - set up of 3100/3100-Avant instruments 27
- emission spectra 17
- equipment, not included in kit 107
- extra peaks, causes 72

F

- fluorescent dyes 16
- FSA sample files 34, 46
- FTA cards
 - amplification 23
 - bloodstained 23

G

GeneMapper™ ID Software
 data analysis 34
overview 16 GeneMapper™
ID-X Software
 data analysis 46
 overview 16
GeneScan size standard
 about 18
 dye label 16
 volume per reaction 28, 29, 31
GS 500 18
GS 600 18

H

hematin, effects of 85
Hi-Di formamide, volume per reaction 28, 29, 31
humic acid, effects of 87

I

identity, probability of 94
import
 HID size standard 42, 55
inheritance 80
instrumentation
 310 genetic analyzer 16, 31
 3100/3100-Avant genetic analyzer 16
 3130/3130xl genetic analyzer 16
 3500/3500xL genetic analyzer 29
 software compatibility 16

K

kit
 allelic ladder 17
 amplification 11
 contents 17
 control DNA 17
 description 11
 fluorescent dyes 16
 loci amplification 12
 master mix 17
 primers 12, 17, 20
 purpose 11
 reagents 17
 supported instruments 11
 thermal cyclers for use with 112
kit performance, comparisons
 DNase I 85
 hematin 85
 humic acid 87

L

Limited Product Warranty 118
LIZ size standard
 about 18
 volume per reaction 28, 29, 31
loci
 characterization 79
 linkage 80
 mapping 80
low TE buffer 19

M

master mix, volume per reaction 21
materials and equipment
 included in kit 17
 not included in kit 107
mixed samples, resolution of genotypes 89
multicomponent analysis 16
mutation studies 93

N

negative control, sample preparation 22
new primers 91

O

off-ladder alleles 64
operating systems 27, 29, 31

P

paternity exclusion, probability of 104
PCR
 components, validation of 60
 cycle number, validation 62
 hematin, inhibitor 85
 humic acid, inhibitor 87
 performing 22
 setup 111
 thermal cycling conditions, programming 22
PCR work areas 111
percent stutter 72
positive control, sample preparation 22
primers
 Amelogenin 79
 volume per reaction 21
probability of identity, definition 103
project examination and editing 58

Q

quantification, DNA 19

R

reaction mix, for PCR 21
reagents, user supplied 19
reproducibility 63
run module, electrophoresis 27, 29, 31

S

safety
 biohazard 114
 chemical 114
sample files, .fsa 34, 46
sample preparation 22
 DNA negative control 22
 DNA positive control 22
 standards 18
sensitivity 82
size deviation, sample alleles and ladder alleles 63
size standard 18
sizing precision 64
software
 overview 45
software, instrument compatibility 16
species specificity 81
split peaks, +A nucleotide addition 76
STRBase 91
stutter 72
stutter percentages, marker-specific 75
support, obtaining 118

T

technical support 118
Terms and Conditions 118
thermal cyclers
 for use with kit 112
thermal cycling
 parameters, validation of 61
 programming conditions 22
training, information on 118

U

user-supplied reagents 19

V

validation
 accuracy, precision, reproducibility 63
 annealing temperatures 61, 62
 characterization of loci 79
 conditions 59
 developmental 60

effect of DNA quantity 83
loci characterization 79
mixture studies 88
mutation rate 93
PCR cycle number 62
population data 91
probability of identity 94
probability of paternity exclusion 104
sensitivity 82
size deviation, sample and ladder alleles 63, 64
species specificity 81
stability 84
stutter 72
thermal cycling parameters 61

W

warranty 118
work area
 amplified DNA 112
 PCR setup 111
 setup and lab design 111
workflow overview 15

