appliedbiosystems

NGM Detect[™] PCR Amplification Kit USER GUIDE

Catalog Number A31832
Publication Number 100044085
Revision D





Life Technologies Ltd | 7 Kingsland Grange | Woolston, Warrington WA1 4SR | United Kingdom For descriptions of symbols on product labels or product documents, go to **thermofisher.com/symbols-definition**.

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. 100044085

Revision	Date	Description
D	2 January 2019	We redesigned primers at the IQC Large and TH01 markers to correct high background seen in samples with a high soil burden, then performed subsequent validation work. Accordingly, we made the following changes to this user guide: Chapter 1—Updated "About the primers", Figure 1, and Figure 2. Chapter 2—Updated Figure 3. Chapter 6—Updated text and figures throughout. Chaper 4—Updated analysis file references to v3.
С	13 March 2017	 Added information for artifact (p 85). Removed IQC from precision and accuracy graphs (update three figures) and tables; noted that they were not included in the study (p 60-66). Updated wording on how stutter filters are calculated; updated minus stutter values; added plus stutter values (pp 75, 81-82). Corrected typo—change 401 to 410 (p 88). Updated SWGDAM references from 2012 to 2016.
В	28 December 2016	 Added Chapter 6, "Experiments and results". Added information about direct amplification to Chapter 2.
А	21 September 2016	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. Windows and Windows Vista are trademarks of Microsoft Corporation. Whatman and FTA are trademarks of Whatman Limited. NUCLEIC-CARD and Copan are trademarks of Copan Italia S.P.A., and used by Life Technologies under their permission. Harris Micro-Punch is a trademark of Harris, Joel S. TA Shunderson Communications. VWR Scientific is a trademark of VWR International, Inc. Robbins Scientific is a trademark of Molecular Bioproducts, Inc. Agilent is a trademark of Agilent Technologies, Inc. Adobe, Acrobat, and Reader are trademarks of Adobe Systems Incorporated.

©2019 Thermo Fisher Scientific Inc. All rights reserved.

Contents

CHAPTER 1 Product information	8
Product description	
Kit overview	
Internal quality control system for PCR	
Validated DNA input amounts and PCR cycles	
About the primers	
Dyes used in the kit	
Loci amplified by the kit	
Standards and controls that are required	
Allelic ladder profile	
DNA Control 007	
Contents and storage	
Required materials not supplied	14
Instruments and software compatibility	15
Workflow	16
CHAPTER 2 Perform PCR	17
DNA quantification	17
Importance of quantification before STR analysis	
Effect of DNA quantity on results	17
Methods of quantifying DNA	18
Before you begin	19
Prepare low-TE buffer	
Thaw reagents (before first use of the kit)	
Prepare the amplification kit reactions	
Perform PCR	
Direct amplification	22

CHAPTER 3 Perform electrophoresis	23
Allelic ladder requirements for electrophoresis	23
Materials required for electrophoresis	24
Set up the 3500/3500xL instruments for electrophoresis (before first use of the kit)	
Electrophoresis software setup	
Create a 3500 QC protocol	
Perform spectral calibration	
Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)	
Set up the 3130/3130 <i>xl</i> instruments for electrophoresis (before first use of the kit)	
Electrophoresis software setup Obtain and activate 6-dye license	
Perform spectral calibration	
Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)	
CHAPTER 4 Analyze data with GeneMapper $^{\text{\tiny TM}}$ <i>ID-X</i> Software	32
Overview of GeneMapper ID-X Software	32
GeneMapper $^{^{ exttt{ iny ID-}X}}$ Software analysis of the NGM Detect $^{^{ exttt{ iny K}}}$ kit	33
Allelic ladder requirements for data analysis	34
File names and versions used in this section	34
Set up the GeneMapper $^{\text{\tiny M}}$ /D-X Software for analysis (before first use of the kit)	35
Workflow: Set up GeneMapper $^{^{\mathrm{TM}}}$ /D-X Software	35
Check panel, bin, and stutter file versions on your computer	
(If needed) Download newer versions of panel, bin, and stutter files	
Import panels, bins, and marker stutter	
(Optional) Define custom table or plot settings	
Create an analysis method	
Create an analysis method	
Enter Analysis Method settings	
Create a size standard definition file if needed	
About the GS600_LIZ_ (60–460) size standard definition file	
Analyze and edit sample files with GeneMapper /D-X Software	
Examine or edit a project	
For more information on using the GeneMapper $^{\text{TD}}$ /D-X Software	
For more information on using the Genemapper ID-X Software	ЭI

CHAPTER 5 Assess the PCR reaction with the Internal Quality Control System	. 52
Overview of the Internal Quality Control System Evaluate the PCR reaction Balanced profile Ski slope profile with decreased IQCL peak height Ski slope profile with balanced IQC peaks No sample peaks with balanced IQC peaks	. 52 . 53 . 54 . 55
CHAPTER 6 Experiments and results	. 56
Importance of validation	. 56
Experiment conditions	. 56
Laboratory requirements for internal validation	57
Developmental validation	. 57
SWGDAM guideline 2.2.1	. 57
SWGDAM guideline 3.9.2	. 57
SWGDAM guideline 3.9.6	. 57
PCR components	. 58
Thermal cycling temperatures	. 58
PCR cycle number	. 60
CE injection time	. 60
Accuracy, precision, and reproducibility	. 63
SWGDAM guideline 3.5	. 63
Accuracy observation	. 63
Precision and size window description	
Precision and size window observation	. 65
Extra peaks in the electropherogram	. 77
Causes of extra peaks	. 77
Extra peaks: Stutter	
Extra peaks: Addition of 3' A nucleotide	. 85
Extra peaks: Artifacts	. 87
Characterization of loci	. 88
SWGDAM guideline 3.1	. 88
Loci in this kit	. 88
Nature of polymorphisms	. 88
Inheritance	. 88
Mapping	. 88
Genetic linkage	. 88
Species specificity	89
SWGDAM Guideline 3.2	
Nonhuman studies	89

Sensitivity
SWGDAM guideline 3.3
Sensitivity observation
Stability 93
SWGDAM guideline 3.4
Degraded DNA
Effect of inhibitors: hematin
Effect of inhibitors: humic acid
Mixture studies 97
SWGDAM guideline 3.8
Mixture study overview
Mixture study observation 98
Resolution of genotypes in mixed samples
Limit of detection of the minor component
Population data
SWGDAM guideline 3.7
Population data overview
Loci in the kit
Population samples used in these studies
Concordance studies
Probability of Identity definition
Probability of identity
Probability of paternity exclusion observation
APPENDIX A Troubleshooting
APPENDIX B Materials required but not supplied
STR kit required materials
Sample preparation required materials
Thermal cycler required materials
ProFlex [™] PCR System
Veriti [™] Thermal Cycler
GeneAmp [™] PCR System 9700
Genetic analyzer required materials
3500 Series Genetic Analyzer
3130 Series Genetic Analyzer required materials
Analysis software required materials
GeneMapper [™] <i>ID-X</i> Software v1.5.2 patch
GeneMapper ID-X Software V1.5.2 patch
Miscellaneous required materials
Plates and tubes
Laboratory supplies

APPENDIX C PCR work areas 1	31
Work area setup and lab design	
PCR setup work area materials	131
Amplified DNA work area	132
APPENDIX D Safety	33
Chemical safety	134
Biological hazard safety	135
Documentation and support	36
Related documentation	136
Customer and technical support	138
Limited product warranty	138
References	39
Index 1	42



Product information

Product description	. 8
Contents and storage	14
Required materials not supplied	14
Instruments and software compatibility	15
Workflow	16

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

Product description

Kit overview

The Applied Biosystems $^{\text{\tiny M}}$ NGM Detect $^{\text{\tiny M}}$ PCR Amplification Kit is a 6-dye, short tandem repeat (STR) multiplex assay for the amplification of specific loci in human genomic DNA.

The kit amplifies:

- 16 autosomal STR loci:
 - The 12 loci in the extended European Standard Set (ESS): FGA, TH01, vWA, D3S1358, D8S1179, D18S51, D21S11, D12S391, D1S1656, D2S441, D10S1248, and D22S1045
 - 3 additional loci that are not in the ESS, but that are present in the AmpFℓSTR[™] SGM Plus[™] PCR Amplification Kit: D16S539, D2S1338, and D19S433
 - SE33, a highly polymorphic locus
- Two internal quality control markers (IQCS and IQCL)
- 1 insertion/deletion polymorphic marker on the Y chromosome (Y indel)
- Amelogenin (sex determining marker)

Internal quality control system for PCR

The NGM DetectTM kit is an Applied BiosystemsTM STR kit that includes an internal quality control (IQC) system for PCR. The IQC system has two synthetic targets, one low molecular weight and one high molecular weight, that are amplified with the sample. The behavior of the IQC target peaks can be used to evaluate the success of the PCR reaction and give an indication of sample quality.

Validated DNA input amounts and PCR cycles

The kit is validated for use with 500 pg DNA (15-µL input volume) for 30 cycles.

About the primers

The NGM Detect^{$^{\text{TM}}$} kit primers are manufactured using the same synthesis and purification improvements as the primers in the GlobalFiler^{$^{\text{TM}}$} kit, Identifiler ^{$^{\text{TM}}$} Plus kit, and NGM SElect^{$^{\text{TM}}$} kit. These improvements enhance the assay signal-to-noise ratio and simplify the interpretation of results.

The primers that are used in the NGM DetectTM kit are unique when compared with primers for corresponding STR loci in the GlobalFilerTM, IdentifilerTM Plus, and NGM SElectTM kits except for one D12 primer and one TH01 primer that are shared with the GlobalFilerTM and NGM SElectTM kits.

Changes to the IQC Large and TH01 primers

A small number of laboratories reported high background with the original formulation of the NGM Detect $^{\text{\tiny{IM}}}$ kit when samples with a high soil burden were run. After internal investigations, we determined that cross-reactivity of the IQC Large and TH01 primers with soil-associated microbial DNA caused the background artifact peaks. Accordingly, we redesigned primers for both markers. We did not make any other changes to the NGM Detect $^{\text{\tiny{IM}}}$ kit formulation, protocols, or workflow.

The updated formulation of the NGM Detect^{$^{\text{M}}$} kit was re-validated with internal and external testing. Other than correcting the soil-associated background issue, the performance of the updated formulation assay is fully comparable to that of the original kit. Unless otherwise indicated, the data in Chapter 6, "Experiments and results" are from the re-validation study.

See the *Technical Note: Updated NGM Detect*[™] *PCR Amplification Kit: Validation and Comparative Study* for studies directly related to soil specificity and a direct comparison of the NGM Detect[™] kit original formulation to the updated formulation.

Dyes used in the kit

Dye	Color	Label
6-FAM [™]	Blue	Samples, allelic ladders, and controls
VIC™	Green	
TED™	Yellow	
TAZ™	Red	
SID [™]	Purple	
LIZ™	Orange	GeneScan [™] -600 LIZ [™] Size Standard v2.0

Chapter 1 Product information Product description

Loci amplified by the kit

Table 1 NGM Detect[™] kit loci and alleles

Locus designation	Chromosome location	Alleles included in Allelic Ladder	Dye label	DNA Control 007
IQCS	N/A	1, 2	6-FAM [™]	2
D2S1338	2q35	11–28		20,23
SE33	6q14	4.2, 6.3, 8, 9, 11–20, 20.2, 21, 21.2, 22.2, 23.2, 24.2, 25.2, 26.2, 27.2, 28.2, 29.2, 30.2, 31.2, 32.2,33.2, 34.2, 35, 35.2, 36–39, 42		17, 25.2
IQCL	N/A	1, 2		2
D16S539	16q24.1	5, 8–15	VIC™	9,10
D18S51	18q21.33	7, 9, 10, 10.2, 11–13, 13.2, 14, 14.2, 15–27		12, 15
TH01	11p15.5	4-9, 9.3, 10-13, 13.3		7, 9.3
D12S391	12p13.2	14–19, 19.3, 20–27		18, 19
D3S1358	3p21.31	9–20	TED™	15, 16
FGA	4q28	13–30, 30.2, 31.2, 32.2, 33.2, 42.2, 43.2, 44.2, 45.2, 46.2, 47.2, 48.2, 50.2, 51.2		24, 26
Y indel	Yq11.221	1, 2	TAZ [™]	2
Amelogenin	X p22.1-22.3, Y: p11.2	X, Y		X, Y
vWA	12p13.31	11–24		14, 16
D21S11	21q11.2-q21	24, 24.2, 25–28, 28.2, 29, 29.2, 30, 30.2, 31, 31.2, 32, 32.2, 33, 33.2, 34, 34.2, 35, 35.2, 36–38		28, 31
D1S1656	1q42.2	9–14, 14.3, 15, 15.3, 16, 16.3, 17, 17.3, 18, 18.3, 19.3, 20.3		13, 16
D2S441	2p14	8–11, 11.3, 12–17		14, 15
D8S1179	8q24.13	5–19	SID [™]	12, 13
D19S433	19q12	6–12, 12.2, 13, 13.2, 14, 14.2, 15, 15.2, 16, 16.2, 17, 17.2, 18.2, 19.2		14, 15
D22S1045	22q12.3	7–20		11, 16
D10S1248	10q26.3	8–19		12, 15

Standards and controls that are required

For the NGM Detect $^{\text{TM}}$ kit, the panel of standards needed for PCR amplification, PCR product sizing, and genotyping are:

- DNA Control 007—A positive control for evaluating the efficiency of the amplification step and STR genotyping using the NGM Detect[™] Allelic Ladder. DNA Control 007 is present in the kit. See "DNA Control 007" on page 13.
- GeneScan[™]-600 LIZ[™] Size Standard v2.0 Used for obtaining sizing results. This standard, which has been evaluated as an internal size standard, yields precise sizing results for PCR products. Order the GeneScan[™]-600 LIZ[™] Size Standard v2.0 (Cat. No. 4408399) separately.
- NGM Detect[™] Allelic Ladder—Developed for accurate characterization of the alleles amplified by the kit. The Allelic Ladder is present in the kit and allows automatic genotyping of most of the reported alleles for the loci in the kit. See "Allelic ladder profile" on page 11.

Allelic ladder profile

The allelic ladder profile appears on the next page.

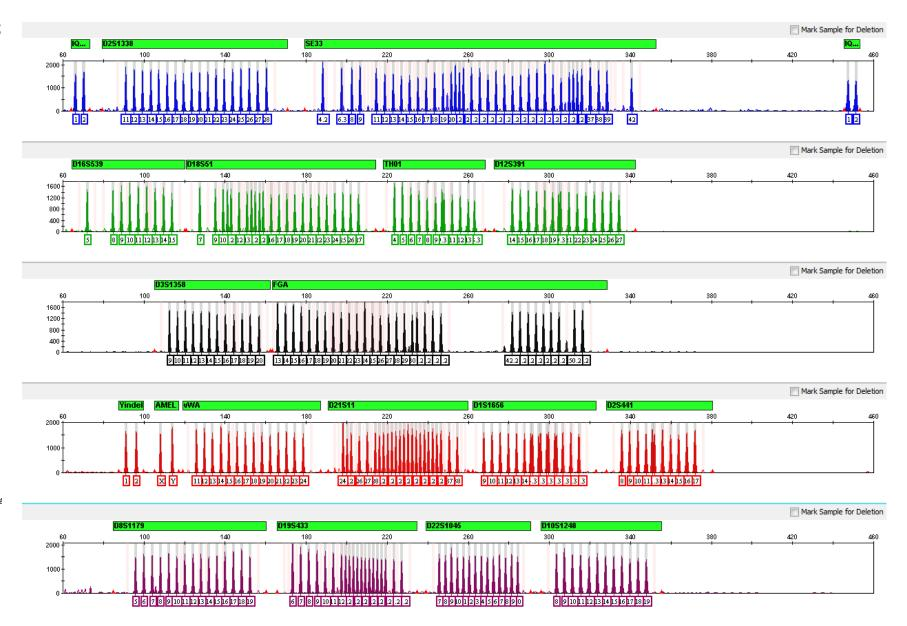


Figure 1 GeneMapper[™] /D-X Software plot of the NGM Detect[™] Allelic Ladder

DNA Control 007

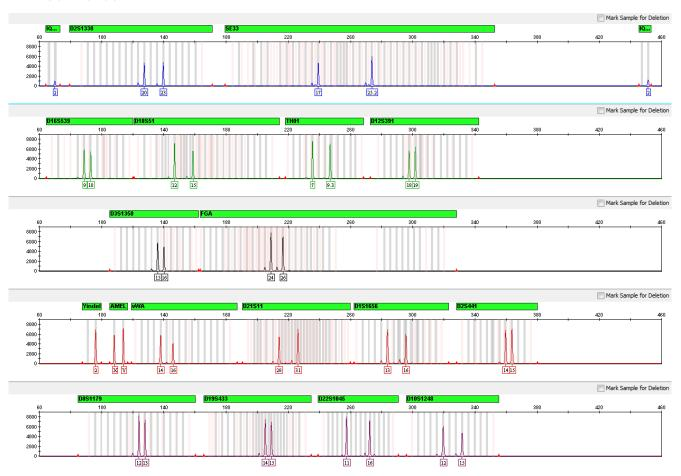


Figure 2 DNA Control 007 (500 pg) amplified with the NGM Detect[™] kit and analyzed on an Applied Biosystems[™] 3500xL Genetic Analyzer (Y-axis scale 0 to 8,000 RFU).

Contents and storage

The NGM Detect $^{\text{\tiny TM}}$ kit (Cat. No. A31832) contains sufficient quantities of the following reagents to perform 200 amplifications with a 25 μ L total reaction volume.

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.

IMPORTANT! Do not refreeze kit components after thawing.

Table 2 NGM Detect[™] PCR Amplification Kit (Cat. No. A31832; 200 reactions)

Contents	Description	Amount	Storage
NGM Detect [™] Master Mix	Contains enzyme, salts, dNTPs, bovine serum albumin, enzyme, and 0.05% sodium azide in buffer and salt.	2 × 0.75 mL	-25°C to -15°C on receipt. 2°C to 8°C after first use, for up to 6 months or up to the expiration date stated on the kit (whichever comes first).
NGM Detect [™] Primer Set	Contains forward and reverse primers to amplify DNA targets.	2 × 0.25 mL	-25°C to -15°C on receipt. 2°C to 8°C after first use, for up to 6 months or up to the expiration date stated on the kit (whichever comes first). Store protected from light.
NGM Detect [™] Allelic Ladder	Contains amplified alleles. See "Allelic ladder profile" on page 11 for information.	1 × 0.05 mL	-25°C to -15°C on receipt. 2°C to 8°C after first use, up to the expiration date stated on the kit. Store protected from light.
DNA Control 007	Contains 0.1 ng/µL human male genomic DNA from cell line in 0.05% sodium azide and buffer ^[1] See "DNA Control 007" on page 13 for information.	1 × 0.3 mL	-25°C to -15°C on receipt. 2°C to 8°C after first use, up to the expiration date stated on the kit.

^[1] DNA Control 007 is included at a concentration that is appropriate for use as an amplification control (that is, to provide confirmation of the capability of the kit reagents to generate a profile of expected genotype). It is not designed for use as a DNA quantification control. If you quantify aliquots of DNA Control 007, the concentration may differ from the labeled concentration.

Required materials not supplied

See Appendix B, "Materials required but not supplied".

Instruments and software compatibility

Instrument type	Validated models		
Thermal	ProFlex [™] 96-well PCR System (Cat. No. 4484075)		
cyclers	ProFlex [™] 2 × 96-well PCR System (Cat. No. 4484076)		
	ProFlex [™] 3 × 32-Well PCR System [Cat. No. 4484073]		
	Veriti [™] 96-Well Thermal Cycler (Cat. No. 4479071)		
	GeneAmp [™] PCR System 9700, 96-Well Silver (Cat. No. N8050001)		
	• GeneAmp [™] PCR System 9700, 96-Well Gold-Plated (Cat. No. 4314878)		
	IMPORTANT! The NGM Detect [™] kit is NOT validated for use with:		
	 ProFlex[™] 2 × Flat PCR System (Cat. No. 4484078) 		
	• ProFlex [™] 2 × 384-well PCR System (Cat. No. 4484077)		
	• Veriti [™] 96-Well <i>Fast</i> Thermal Cycler (Cat. No. 4375305)		
	• GeneAmp [™] PCR System 9700 with the aluminium 96-well block (Cat. No. 4314879)		
Genetic	3500/3500xL Genetic Analyzer with any of the following:		
analyzers ^[1]	 3500 Data Collection Software v1 (Windows[™] Vista operating system) and HID Updater 3500 Data Collection Software v2 (Cat. No. 4480670) 		
	 3500 Data Collection v2 Software (Windows[™] 7 operating system) and HID Updater 3500 Data Collection Software v2 (Cat. No. 4480670) 		
	 3500 Data Collection v3 Software (Windows[™] 7 operating system) 		
	• 3130/3130 <i>xl</i> Genetic Analyzer with:		
	 Data Collection Software v4 (Windows[™] 7 operating system) 		
	- 3130/3730 Data Collection v4 6-Dye Module v1		
Analysis	GeneMapper [™] ID - X Software v1.5.2 or later ^[2]		
software	Windows [™] 7 operating system		

 $^{^{[1]}}$ We conducted validation studies using the 3130xl, 3500, and 3500xL configurations.

^[2] GeneMapper[™] /D-XSoftware v1.2 to v1.5 can be used to analyze NGM Detect[™] PCR Amplification Kit data. However, some genotype quality assessment features of the NGM Detect[™] kit are not included in earlier versions of the software. Refer to Chapter 4, "Analyze data with GeneMapper[™] ID-X Software" for more details.

Workflow

Extract DNA, see:

www.thermofisher.com/hid-sampleprep



Quantify DNA

"DNA quantification" on page 17



Perform PCR

"Prepare the amplification kit reactions" on page 19



"Perform PCR" on page 21



Perform electrophoresis

"Set up the 3500/3500xL instruments for electrophoresis (before first use of the kit)" on page 24 or "Set up the 3130/3130xl instruments for electrophoresis (before first use of the kit)" on page 28



"Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)" on page 27



Analyze data

"Set up the GeneMapper $^{\text{\tiny M}}$ ID-X Software for analysis (before first use of the kit)" on page 35



"Create an analysis method" on page 40



"Create a size standard definition file if needed" on page 48



"Analyze and edit sample files with GeneMapper ID-X Software" on page 50



"Examine or edit a project" on page 51



Chapter 5, "Assess the PCR reaction with the Internal Quality Control System"



Perform PCR

DNA quantification	17
Before you begin	19
Prepare the amplification kit reactions	19
Perform PCR	21
Direct amplification	22

DNA quantification

Importance of quantification before STR analysis

DNA quantification can be used to determine:

- If the sample contains sufficient human DNA and/or human male DNA to proceed with short tandem repeat (STR) analysis.
- The amount of sample to use in STR analysis applications.
- The relative quantities of human male and female DNA in a sample (using the Quantifiler[™] Trio DNA Quantification Kit). This guides selection of the applicable STR chemistry.
- The DNA quality, with respect to the inhibition level and the DNA degradation level. This metric is useful for determining the likelihood of recovery of STR loci with larger amplicon sizes.
- If the sample contains highly degraded DNA. Such samples may require an
 alternative approach to STR analysis by capillary electrophoresis. Precision ID
 NGS System and Panels are optimized for degraded samples. The Precision ID
 Identity Panel provides discrimination of individuals similar to STR genotype
 match probabilities. Also, the Precision ID Ancestry Panel infers biogeographical
 ancestry for investigative leads.
- If PCR inhibitors are present in a sample. Such that samples may require additional purification before proceeding to STR analysis.

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 capillary electrophoresis instrument ("off-scale" data).

Chapter 2 Perform PCR DNA quantification

Off-scale data are 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 a corresponding stutter peak to appear higher in relative intensity, therefore increasing the calculated percent stutter.
- Multicomponent analysis of off-scale data are not accurate. This inaccuracy results in poor spectral separation ("pull-up").
- False signals of inhibition by the IQC system, although none is present.
- A reduction in the IQCL peak height.
- Incomplete +A nucleotide addition.

To address these problems, rerun the amplification reaction using less DNA.

If too little DNA is added to the PCR reaction, the total number of allele copies added to the PCR is extremely low. Unbalanced amplification of the alleles may occur because of stochastic fluctuation.

Methods of quantifying DNA

For information on recent innovations in quantification chemistry, go to **thermofisher.com**.

Kit and user guide	Detects	How it works
Quantifiler [™] HP DNA Quantification Kit (Cat. No. 4482911) For more information, see Quantifiler [™] HP and Quantifiler [™] Trio DNA Quantification Kits User Guide (Pub. No. 4485354)	 Total human DNA (two targets—one small amplicon and one larger amplicon) Degraded DNA 	 Uses 5' nuclease assays with multiple-copy target loci, for improved detection sensitivity:^[1] The human-specific target loci are multiple copy, and dispersed on various autosomal chromosomes. The primary quantification targets have relatively short amplicons (75 to 80 bases), to improve the detection of degraded DNA samples.
Quantifiler [™] Trio DNA Quantification Kit (Cat. No. 4482910) For more information, see Quantifiler [™] HP and Quantifiler Trio DNA Quantification Kits User Guide (Pub. No. 4485354)	 Total human DNA (two targets—one small amplicon and one larger amplicon) Human male DNA Degraded DNA 	 Uses features that maximize consistency of quantification: Genomic targets have conserved primer- and probe-binding sites. Minimal copy number variation between different individuals and population groups. Contains a Large Autosomal target with a longer amplicon (>200 bases) to help determine if a DNA sample is degraded. Contains an Internal PCR control (IPC) 5' nuclease assay which amplifies an integrated synthetic DNA sequence. The performance of this assay can be used to assess whether real-time PCR of the sample has been impacted by inhibition.

^[1] The detection sensitivity of the Quantifiler™ HP Kit and the Quantifiler™ Trio Kit is improved over the Quantifiler™ Duo Kit.

Before you begin

Prepare low-TE buffer

For optimal results, we recommend using low-TE buffer for sample preparation. Prepare it as described in this procedure or buy it from Teknova (Cat. No. T0223).

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

Note: Adjust the volumes accordingly for specific needs.

- 2. Aliquot, then autoclave the solutions.
- **3.** Store the aliquots at room temperature.

Thaw reagents (before first use of the kit)

Thaw the Master Mix and Primer Set.

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.

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

Prepare the amplification kit reactions

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.

IMPORTANT! Use adhesive film for plate sealing to provide a consistent seal across all wells and prevent evaporation. Caps may not provide a consistent seal across all plate wells.

- 1. Vortex the Master Mix and Primer Set for 3 seconds. Before opening the tubes, remove droplets from the caps by centrifuging the tubes briefly.
- 2. Pipette the required volumes of components into an appropriately sized clear (non-colored) polypropylene tube:

Reaction component	Volume per reaction
Master Mix	7.5 μL
Primer Set	2.5 μL

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

- **3.** Vortex the reaction mixture for 3 seconds, then centrifuge briefly.
- **4.** Dispense 10 μ L of reaction mixture into each reaction well of a MicroAmpTM Optical 96-Well Reaction Plate or each MicroAmpTM tube.
- **5.** Adjust the sample input amount and volume as needed:

Note: The optimum DNA input is 500 pg.

- If total sample input amount is >500 pg of DNA, dilute with low-TE buffer to achieve a 500 pg input in a 15-μL volume.
- If total sample input volume is <15 μ L, bring to volume with low-TE buffer to achieve a 15- μ L input volume.
- **6.** Prepare samples as shown in the following table, then add them to the appropriate well or tube (final reaction volume is $25 \,\mu L$).

Sample	Add		
Negative control	15 μL of low-TE buffer		
Test sample	15 μL of DNA ^[1]		
Positive control	Combine, then add to the reaction well or tube: • 5 µL DNA Control 007 (0.1 ng/µL) • 10 µL of low-TE buffer		

^[1] Prepared in step 5.

- 7. Seal the MicroAmp[™] Optical 96-Well Reaction Plate with MicroAmp[™] Clear Adhesive Film or MicroAmp[™] Optical Adhesive Film.
- **8.** Vortex the plate or tubes at medium speed for 3 seconds.
- **9.** Centrifuge the tubes or plate at 3,000 rpm for approximately 30 seconds in a tabletop centrifuge (with plate holders, if using 96-well plates).
- **10.** Amplify the samples.

IMPORTANT! See "Instruments and software compatibility" on page 15 for a list of validated thermal cyclers.

Perform PCR

IMPORTANT! This kit is validated for use with the thermal cyclers listed in "Instruments and software compatibility" on page 15.

1. Program the thermal cycling conditions.

IMPORTANT! If you are using the:

- ProFlex[™] PCR System, select 9700 Simulation Mode.
- GeneAmp[™] PCR System 9700, select the Max ramping mode.
- Veriti[™] Thermal Cycler, set up the method using the Convert a Method tool and select 9700 Max Mode.

Do not use 9600 emulation mode.

Initial	Cycle (30 cycles)			Final	-
incubation step	Denature Anneal Exte		Extend	extension	Final hold
HOLD	CYCLE			HOLD	HOLD
95°C, 1 minute	96°C, 59°C, 5 seconds 21 seconds		65°C, 29 seconds	60°C, 5 minutes	4°C ^[1]

^[1] The infinity (∞) setting allows an unlimited hold time.

2. Load the plate into the thermal cycler, close the heated cover, then start the run.

IMPORTANT! If you are using a GeneAmp[™] PCR System 9700 *and* adhesive clear film instead of caps to seal the plate wells, place a MicroAmp[™] Optical Film Compression Pad (Cat. No. 4312639) on top of the plate to prevent evaporation during thermal cycling. The Veriti[™] Thermal Cycler and the ProFlex[™] PCR System do not require a compression pad.

3. When the run is complete, store the amplified DNA.

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

IMPORTANT! Protect the amplified DNA from light.

Direct amplification

 $\mathsf{FTA}^{\mathsf{TM}}$ cards and non-chemically treated bloodstain cards are 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 or 96-well plate, washed, then amplified, without transferring the disc.

Our studies indicate that a 1.2-mm bloodstained disc contains approximately 5 ng—20 ng of DNA. Because of the high quantity of DNA, a lower cycle number is required to produce on-scale data. In our testing, an appropriate cycle number for this high quantity DNA was 26 cycles. We recommend that each laboratory determine the optimum cycle number which is based on internal validation studies.

Note: This kit is not fully validated for use with direct amplification. Perform your own validation for this purpose or use the NGM $SElect^{TM}$ Express kit for direct amplification of database samples.

In the example that is shown in Figure 3, a 1.2-mm disc of a bloodstained FTATM card was purified using one wash with FTATM Purification Reagent and one wash with 1X low TE buffer, followed by a short drying step. The sample was then amplified directly in the well of a standard 96-well amplification MicroAmpTM plate for 26 cycles.

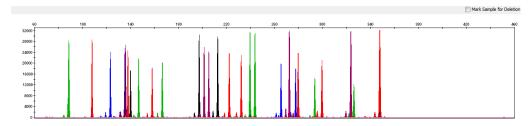


Figure 3 Combined dyes electropherogram of a 1.2-mm FTA[™] bloodstain disc amplified for 26 cycles with the NGM Detect[™] kit on an Applied Biosystems[™] 3500xL Genetic Analyzer [Y-axis scale 0 to 33,000 RFU].



Perform electrophoresis

Allelic ladder requirements for electrophoresis	23
Materials required for electrophoresis	24
Set up the 3500/3500xL instruments for electrophoresis (before first use of the kit)	24
Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)	27
Set up the 3130/3130xl instruments for electrophoresis (before first use of the kit)	28
Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)	30

Allelic ladder requirements for electrophoresis

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

Instrument	Number of allelic ladders to run	One injection equals	Number of samples per allelic ladder(s)
3500	1 per 3 injections	8 samples	23 samples + 1 allelic ladder
3500 <i>xl</i>	1 per injection	24 samples	23 samples + 1 allelic ladder
3130	1 per 4 injections	4 samples	15 samples + 1 allelic ladder
3130 <i>xl</i>	1 per injection	16 samples	15 samples + 1 allelic ladder

IMPORTANT! Variation in laboratory temperature can cause changes in fragment migration speed and sizing variation between runs. Follow the guidelines in the preceding table, which should account for normal variation in run speed. Perform internal validation studies to 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. Size values obtained for the same sample can differ between instrument platforms, because of different polymer matrices and electrophoretic conditions.

Materials required for electrophoresis

Appendix B, "Materials required but not supplied" lists the required materials that are not supplied with this 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.

Set up the 3500/3500xL instruments for electrophoresis (before first use of the kit)

Electrophoresis software setup

The following table lists the data collection software and the run modules that you can use to analyze PCR products that are generated by this kit. For details on the procedures, see the documents that are listed in "Documentation and support" on page 136.

Note: We conducted validation studies for the kit using the 3130*xl*, 3500, or 3500*x*L configurations.

Genetic Analyzer	Operating System	Data Collection Software	Additional software	Instrument protocols, run modules, and conditions
3500 3500xL	Windows [™] Vista	3500 Data Collection Software v1	HID Updater 3500 DC v2 (Cat. No. 4480670)	 Set up the following conditions: Run module: HID36_P0P4 (HID36_P0P4xl for 3500xL) Injection conditions^[1]: 1.2 kV/11 sec (20 sec for 3500xL) Run conditions: 13 kV/1500 sec (13 kV/1500 sec for 3500xL) Dye Set J6-T
3500 3500xL	Windows [™] 7	3500 Data Collection Software v2	HID Updater 3500 DC v2 (Cat. No. 4480670)	Same as 3500 Data Collection Software v1 listed above
3500 3500xL	Windows [™] 7	3500 Data Collection Software v3	None	Same as 3500 Data Collection Software v1 listed above

Our studies indicate that the injection conditions that are documented generate profiles from 0.5 ng of input DNA with heterozygous peak height averages between 4,000–10,000 RFU (3500 or 3500xL) with no instances of allelic dropout and minimal occurrence of off-scale allele peaks. However, individual CE instrument signal intensities can vary, therefore changes to injection parameters may need to be explored and validated to deliver the best results on your specific system. Large deviations from the recommended injection parameters could affect the performance of the size standard and/or allelic ladder, therefore validation is recommended.

Create a 3500 QC protocol

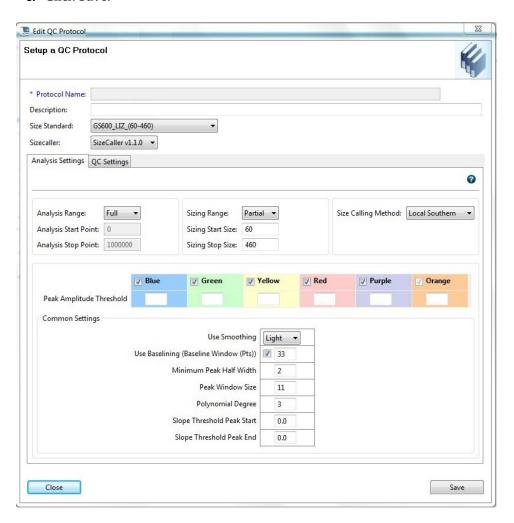
The NGM Detect $^{\text{\tiny TM}}$ kit has been validated with data that was analyzed using the Local Southern method (60–460 base pairs).

- 1. In the Library tab, open the QC Protocol window.
- **2.** Create a new QC protocol:
 - **a.** Name the new QC protocol according to your laboratory naming convention.
 - **b**. Set the following parameters:

Parameter	Setting
Size Standard	GS600_LIZ_(60-460)
Size Range	Partial
Sizing Start Size	60 bp
Sizing Stop Size	460 bp
Size Calling Method	Local Southern Method
After checking the "Use Baselining" box: Baseline Window Pts.	33
Peak Window Size	11



c. Click Save.



3. Add the QC protocol to the HID assay.

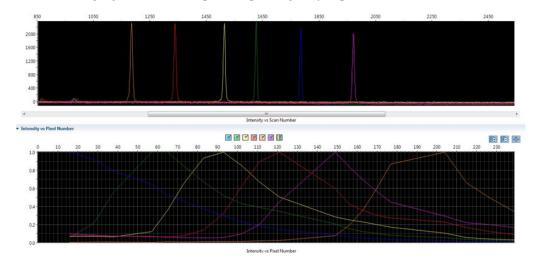
Perform spectral calibration

Perform a spectral calibration using the DS-37 Matrix Standard Kit (Dye set J6-T, 6-dye) (J6-T Dye Set) (Cat. No. A31234).

Note: Since J6-T is a new dye set, create a new dye set in the Data Collection Software before running a spectral calibration. For instructions on creating a new dye set, see the "Create a New Dye Set" section of the 3500/3500xL Genetic Analyzer with 3500 Series

Data Collection Software v2 User Guide (Cat. No. 4476988). Use the J6 template to set up the J6-T dye set.

The following figure is an example of a passing 6-dye spectral calibration.



Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)

This procedure applies to the 3500 Series and 3130 Series instruments.

Prepare the samples for electrophoresis immediately before loading.

1. Pipet the required volumes of components into an appropriately sized polypropylene tube:

Reagent	Volume per reaction
GeneScan [™] -600 LIZ [™] Size Standard v2.0	0.4 μL
Hi-Di [™] Formamide	9.6 µL

Note: Include volume for additional samples 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.** Vortex the tube, then briefly centrifuge.
- 3. 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 10 μL of Hi-Di[™] Formamide.

- 4. 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.
- **5**. Heat the reaction plate in a thermal cycler at 95°C for 3 minutes.
- **6.** Immediately place the plate on ice for 3 minutes.
- 7. Place the sample tray on the autosampler, then start the electrophoresis run.

Set up the 3130/3130*xl* instruments for electrophoresis (before first use of the kit)

Electrophoresis software setup

The following table lists the data collection software and the run modules that you can use to analyze PCR products generated by this kit. For details on the procedures, see the documents listed in "Documentation and support" on page 136.

Genetic Analyzer	Operating System	Data Collection Software	Additional software	Run modules and conditions
3130 ^[1]	Windows [™] 7	Data Collection Software v4	3130/3730 DC v4 6- Dye Module v1	Set up the following conditions: • HIDFragmentAnalysis36_P0P4_1 • Injection conditions ^[2] : 3 kV/4 sec • Run conditions: 14 kV/1600 sec • Dye Set J6-T
3130 <i>xl</i>				Set up the following conditions: • HIDFragmentAnalysis36_P0P4_1 • Injection conditions ^[2] : 3 kV/5 sec • Run conditions: 14 kV/1600 sec • Dye Set J6-T

 $^{^{[1]}}$ We conducted validation studies using the 3130xL, 3500, and 3500xL configurations.

Obtain and activate 6-dye license

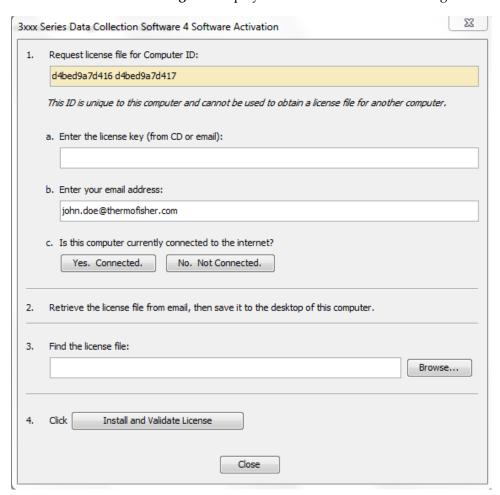
- 1. Confirm that you are running Data Collection Software v4 (Help > About).
- **2.** Obtain a 3130 DC v4 6-Dye Module v1 License key. Contact your local Human Identification representative for information.

^[2] Our studies indicate that the injection conditions that are documented generate profiles from 0.5 ng of input DNA with heterozygous peak height averages between 2,000–4,000 RFU (3130 or 3130xl) with no instances of allelic dropout and minimal occurrence of off-scale allele peaks. However, individual CE instrument signal intensities can vary, therefore changes to injection parameters may need to be explored and validated to deliver the best results on your specific system. Large deviations from the recommended injection parameters could affect the performance of the size standard and/or allelic ladder, therefore validation is recommended.

3. Ensure that all network cards in the computer are enabled.

IMPORTANT! You can run the 3130 Series Data Collection Software v4 using only the network cards that are enabled when you activate the software license. For example, if you activate the software when your wireless network card is disabled, you will not be able to run the software when the wireless network card is enabled.

4. Select Tools > License Manager to display the Software Activation dialog box.



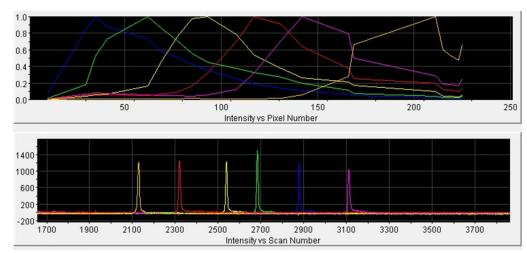
- **5.** Request the software license file by performing steps 1a, 1b, and 1c as listed on the activation screen. The license file will be emailed to you.
- **6.** Obtain the software license file from your email.
- 7. Make a copy of the software license file and keep it in a safe location.
- **8.** Copy the software license file to the desktop of the Data Collection Software v4 computer.
- **9.** If the Software Activation dialog box has closed, select **Tools > License Manager**.

- **10.** Click **Browse**, then navigate to the software license file saved on your computer.
- Click Install and Validate License.
 A message is displayed when the license is installed and validated.
- 12. Click Close.

Perform spectral calibration

Perform a spectral calibration using the DS-37 Matrix Standard Kit (Dye set J6-T, 6-dye) (J6-T Dye Set) (Cat. No. A31234).

The following figure is an example of a passing 6-dye spectral calibration.



Prepare samples for electrophoresis (3500 Series and 3130 Series instruments)

This procedure applies to the 3500 Series and 3130 Series instruments.

Prepare the samples for electrophoresis immediately before loading.

1. Pipet the required volumes of components into an appropriately sized polypropylene tube:

Reagent	Volume per reaction
GeneScan [™] -600 LIZ [™] Size Standard v2.0	0.4 μL
Hi-Di [™] Formamide	9.6 μL

Note: Include volume for additional samples 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. Vortex the tube, then briefly centrifuge.

- **3.** 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 10 μ L of Hi-DiTM Formamide.

- **4.** 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.
- **5.** Heat the reaction plate in a thermal cycler at 95°C for 3 minutes.
- **6.** Immediately place the plate on ice for 3 minutes.
- **7.** Place the sample tray on the autosampler, then start the electrophoresis run.



Analyze data with GeneMapper[™] *ID-X* Software

Overview of GeneMapper ID-X Software	32
$GeneMapper^{^{TM}}ID\text{-X}Software\;analysis\;of\;the\;NGM\;Detect^{^{TM}}kit\;\ldots\ldots$	33
Allelic ladder requirements for data analysis	34
File names and versions used in this section	34
Set up the GeneMapper [™] ID-X Software for analysis (before first use of the kit)	35
Create an analysis method	40
Create a size standard definition file if needed	48
Analyze and edit sample files with GeneMapper ID-X Software	50
Examine or edit a project	51
For more information on using the GeneMapper [™] ID-X Software	51

Overview of GeneMapper ID-X Software

GeneMapper $^{\text{TM}}$ *ID-X* Software is an automated genotyping software application for forensic casework, databasing, and paternity data analysis.

GeneMapperTM ID-X Software v1.5.2 or later analyzes 4-dye, 5-dye, and 6-dye data and is required to correctly analyze data that is generated using the NGM DetectTM kit. After electrophoresis, the data collection software stores information for each sample in a .fsa or .hid file. The GeneMapperTM ID-X Software v1.5.2 or later allows you to analyze and interpret the data from the .fsa or .hid files.

Prior versions of GeneMapperTM ID-X Software (v1.2 and above) are capable of analyzing data that is generated using the NGM DetectTM kit. However, some data quality assessment tools are not fully functional on versions prior to v1.5.2. Details on the data assessment functionality added to specific versions of GeneMapperTM ID-X Software are found in the release notes for each version. To summarize:

- In versions prior to v1.4, samples show a red flag for the CGQ if the Y-indel peak is missing, for example, with a female DNA sample. This occurs because the absence of the Y-indel peak triggers the Allele Number Rule.
- In versions prior to v1.5.2, when the sample type is set as "Negative Control" the software will assign red CGQ flags. This occurs even if the negative control sample is valid, that is, it contains no peaks above the PAT in the STR, AMEL, and Y-indel marker ranges. The red flags appear because the presence of the IQC peaks triggers the Allele Number Rule for this sample type (no peaks are expected in a valid negative control sample).

4

GeneMapper[™] *ID-X* Software analysis of the NGM Detect[™] kit

GeneMapper[™] *ID-X* Software allows you to designate four sample types:

- Sample
- Positive Control
- Allelic Ladder
- Negative Control

GeneMapper $^{\text{TM}}$ *ID-X* Software has logic for Negative Control sample type assignments, whereby the software generates red flags on the Genotype Quality (GQ) indicators when peaks are present above the Peak Amplitude Threshold (PAT) designated in the analysis method.

The NGM Detect[™] PCR Amplification Kit includes the internal quality control (IQC) system for PCR. In this system, two synthetic targets in the primer mix are amplified with the sample under test. The IQC system allows you to determine whether PCR has been successful in the absence of peaks from the sample.

In versions prior to v1.5.2, when the sample type is set as "Negative Control" the software will assign red CGQ flags. This occurs even if the negative control sample is valid, that is, it contains no peaks above the PAT in the STR, AMEL, and Y-indel marker ranges. The red flags appear because the presence of the IQC peaks triggers the Allele Number Rule for this sample type (no peaks are expected in a valid negative control sample).

Thermo Fisher Scientific has developed the NGM Detect^M kit using version 1.5.2 of the GeneMapper^M ID-X Software. This version includes updated logic for the Negative Control sample type. It allows the IQC marker GQ to flag green in the presence of the IQC peaks.

A software patch that updates GeneMapper^{$^{\text{TM}}$} *ID-X* Software v1.5 to v1.5.2 is available at thermofisher.com/us/en/home/technical-resources/software-downloads/genemapper-id-x-software.html.

Allelic ladder requirements for data analysis

- HID analysis requires at least one allelic ladder sample per run folder. Perform
 the appropriate internal validation studies before you use multiple allelic ladder
 samples in an analysis.
 - For multiple allelic ladder samples, the GeneMapper^{$^{\text{TM}}$} *ID-X* Software calculates allelic bin offsets by using an average of all allelic ladders that use the same panel in 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 ladders in 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. Analysis will fail if the Allelic Ladder Sample Type is not specified.
- 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 allelic ladders do exist. Off-ladder (OL) alleles can
 contain full and/or partial repeat units. An off-ladder allele is an allele that occurs
 outside the 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.

File names and versions used in this section

The file names and version numbers of panel, bin, and stutter files that are shown in this section may differ from the file names that you see when you download or import files.

If you need help to determine the correct files to use, contact your local Human Identification representative, or go to **thermofisher.com/support**.

Set up the GeneMapper $^{\text{\tiny TM}}$ /D-X Software for analysis (before first use of the kit)

Workflow: Set up GeneMapper[™] /D-XSoftware Before you use GeneMapper^{$^{\text{TM}}$} *ID-X* Software to analyze data for the first time, you must do the following:

"Check panel, bin, and stutter file versions on your computer" on page 35



"(If needed) Download newer versions of panel, bin, and stutter files" on page 35



"Import panels, bins, and marker stutter" on page 36



"(Optional) Define custom table or plot settings" on page 39

Check panel, bin, and stutter file versions on your computer

- 1. Start the GeneMapper $^{\text{\tiny TM}}$ *ID-X* Software , then log in with the appropriate user name and password.
- 2. Select Tools ▶ Panel Manager.
- 3. Check the version of files that are currently available in the **Panel Manager**:
 - a. Select **Panel Manager** in the navigation pane.
 - **b.** Expand the **Panel Manager folder** and any subfolders to identify the analysis file version that is already installed for your kit choice.
- 4. Check the version of files available for import into the **Panel Manager**:
 - a. Select Panel Manager, then select File ▶ Import Panels to open the Import Panels dialog box.
 - **b.** Navigate to, then open the **Panels** folder, then check the version of panel, bin, and stutter files installed.
- **5.** Check for newer versions of the files as described in the next procedure.

(If needed)
Download newer
versions of panel,
bin, and stutter
files

- 1. Go to thermofisher.com/us/en/home/technical-resources/software-downloads/genemapper-id-x-software.html.
- **2.** If the file versions listed are newer than the versions on your computer, download the file **NGM Detect Analysis Files**.

Note: When downloading new versions of analysis files, see the associated **Read Me** file for details of changes between software file versions. Perform the appropriate internal validation studies before using new file versions for analysis.

3. Unzip the file.



Import panels, bins, and marker stutter

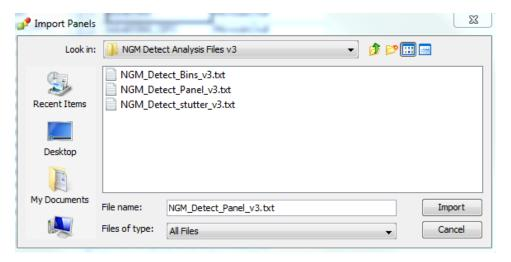
To import the latest panel, bin set, and marker stutter from the website into the GeneMapper $^{\text{TM}}$ *ID-X* Software database:

- 1. Start the GeneMapper^{$^{\text{TM}}$} *ID-X* Software, then log in with the appropriate user name and password.
- 2. Select Tools ▶ Panel Manager.
- **3.** Find, then open the folder containing the panels, bins, and marker stutter:



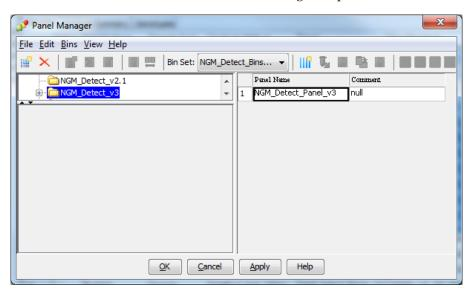
- a. Select Panel Manager, then select File > Import Panels to open the Import Panels dialog box.
- **b.** Navigate to, then open the **NGM Detect Analysis Files** folder that you unzipped in the previous procedure.
- 4. Select NGM_Detect_Panels.txt, then click Import.

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



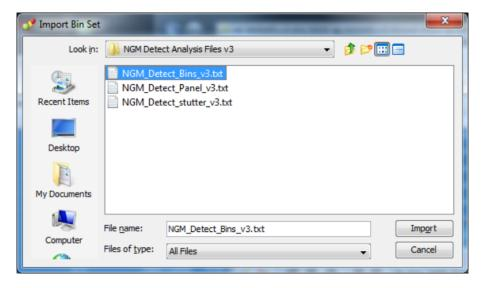
5. Import the bins file:

a. Select the NGM_Detect_Panel 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 Detect Analysis Files folder.
- d. Select NGM_Detect_Bins.txt, then click Import.

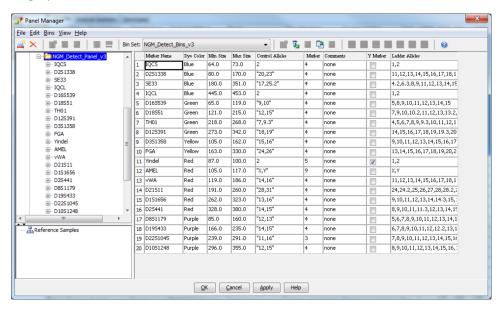
Note: Importing this file associates the bin set with the panels in the **NGM_Detect_Panel** folder.





6. (*Optional*) View the imported panels and bins in the navigation pane: Double-click the **NGM_Detect_Panel** folder.

The panel information is displayed in the right pane and the markers are displayed below it.

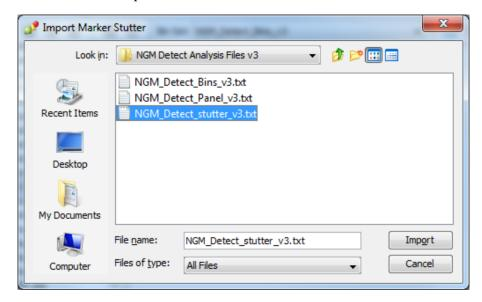


- **7.** Import the stutter file:
 - a. Select the NGM_Detect_Panel folder in the navigation panel.
 - b. Select File > Import Marker Stutter to open the Import Marker Stutter dialog box.
 - c. Navigate to, then open the NGM Detect Analysis Files folder.

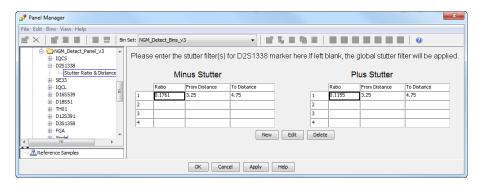
38

d. Select NGM_Detect_Stutter.txt, then click Import.

Note: Importing this file associates the marker stutter ratio with the bin set in the **NGM_Detect_Panel** folder and overwrites any existing stutter ratios associated with the panels and bins in that folder.



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



9. Click **Apply**, then click **OK** to add the panel, bin set, and marker stutter to the GeneMapper[™] *ID-X* Software database.

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

(Optional) Define custom table or plot settings

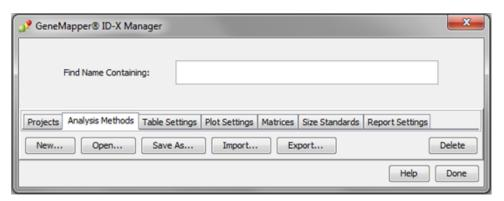
Default views for table and plot settings are provided with the software. For information on defining custom views, see $GeneMapper^{TM}$ ID-X Software Getting Started Guide — Basic Features.

Create an analysis method

Create an analysis method

IMPORTANT! Analysis methods are version-specific, so you must create an analysis method for each version of the software. For example, an analysis method that is created in GeneMapper $^{\text{TM}}$ *ID-X* Software version 1.2 is not compatible with analysis methods that are created in earlier versions of software, or with GeneMapper Software v3.2.1.

 Select Tools ➤ GeneMapper® ID-X Manager to open the GeneMapper ID-X Manager.



- 2. Click 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.

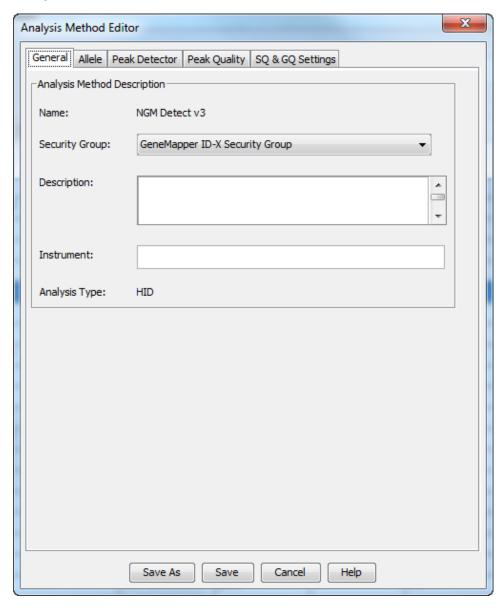
Note: The **Analysis Method Editor** closes when you save your settings. To complete this step quickly, do not save the analysis method until you finish entering settings in all of the tabs.

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

Enter Analysis Method settings

Enter General tab settings

1. Enter a **Name** and select the **Security Group** appropriate for your software configuration.



2. (Optional) Enter a Description and Instrument.

Enter Allele tab settings

IMPORTANT! Perform appropriate internal validation studies to determine the appropriate settings to use.

1. Select the NGM_Detect_Bins_v3 bin set.

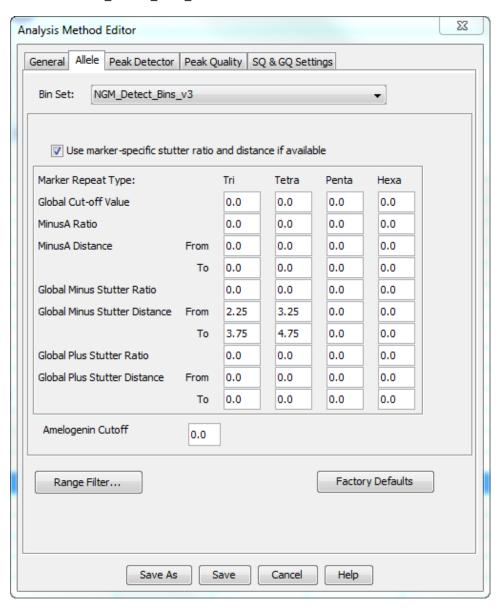


Figure 4 Settings used in developmental validation of the kit

(Optional) To apply the stutter ratios contained in the NGM_Detect_Stutter.txt, select the Use marker-specific stutter ratio and distance if available checkbox (selected by default).

- **3.** If using GeneMapper $^{\text{\tiny TM}}$ *ID-X* Software v1.0.1 or later, enter values for the 4 Marker Repeat Types.
- **4.** Enter the appropriate filter settings.

Enter Peak Detector tab settings

Enter the appropriate values:

Field	Values to enter or select	Additional information
Ranges	Enter the values shown in Figure 5.	_
	Note: The read region for the NGM Detect [™] kit is 64 to 458 bp.	
Peak Detection	Enter the appropriate settings. IMPORTANT! Perform appropriate internal validation studies to determine the appropriate peak amplitude thresholds for interpretation of data.	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.

Field	Values to enter or select	Additional information
Smoothing and Baseline	Enter the values shown in Figure 5, or adjust as needed dependent on the polymer you are using.	_
Size Calling Method	Select Local Southern Method , or another method if validated by your internal validation.	_
Normalization	(Optional) Select the Normalization checkbox.	A Normalization checkbox is available on this tab in GeneMapper [™] / <i>D-X</i> Software for use in conjunction with data run on the 3500 Series Genetic Analyzers.

23 Analysis Method Editor General Allele Peak Detector Peak Quality SQ & GQ Settings Peak Detection Algorithm: Advanced Ranges Peak Detection Analysis Sizing Peak Amplitude Thresholds: Full Range Partial Sizes R: Start Size: 60 Start Pt: 0 G: P: Stop Pt: 10000 Stop Size: 460 Y: 0: Smoothing and Baselining 2 pts Min. Peak Half Width: Smoothing None Polynomial Degree: 3 Light Peak Window Size: 11 pts Heavy Slope Threshold Baseline Window: 33 pts 0.0 Peak Start: Size Calling Method Peak End: 0.0 2nd Order Least Squares 3rd Order Least Squares Normalization Cubic Spline Interpolation Use Normalization, if applicable Local Southern Method Global Southern Method Factory Defaults Save As Save Cancel Help

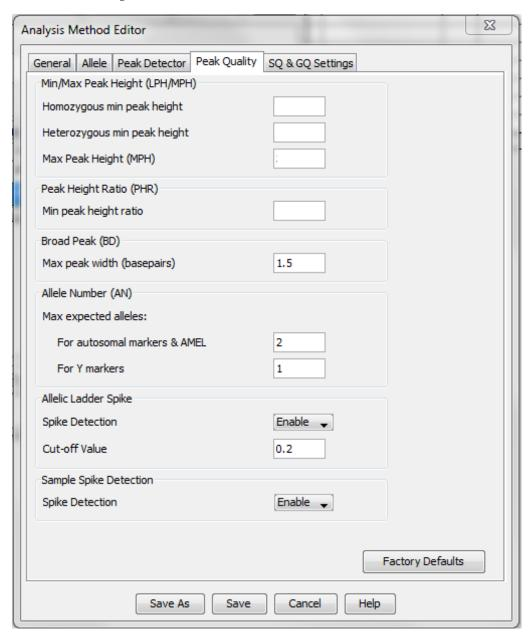
Note: The read region for the NGM Detect[™] kit is 64 to 458 bp.

Figure 5 Settings used in developmental validation of the kit

Enter Peak Quality tab 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 data.

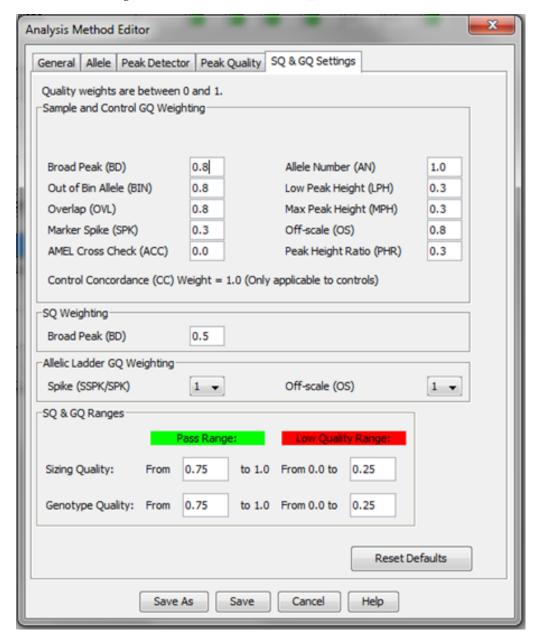
Enter the following values:



Enter SQ and GQ tab settings

IMPORTANT! The values that are 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.

Enter the following values:



Note: Set the **ACC GQ Weighting** according to the values you determine during internal validation studies of the **ACC PQV**. For example, set the **ACC GQ Weighting** to 0.3 or higher to flag samples in which the Amelogenin result is anything other than X, X or X, Y, or does not agree with the results for the Y indel marker.

Create a size standard definition file if needed

If you cannot use the default settings that are provided, create a new size standard definition file.

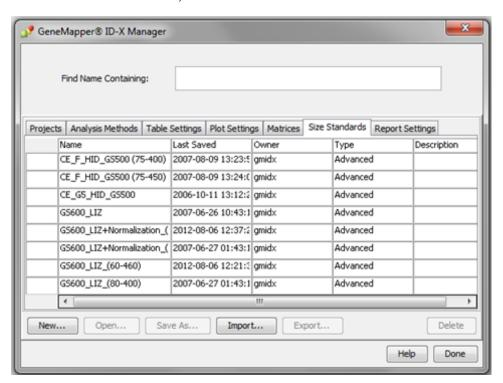
About the GS600_LIZ_ (60-460) size standard definition file

The GS600_LIZ_(60–460) size standard definition that is provided with GeneMapper TD - T

This size standard definition has been validated for use with this kit on the genetic analyzers listed in "Instruments and software compatibility" on page 15. If you need to create your own size standard definition, see "Create a size standard definition file" on page 48.

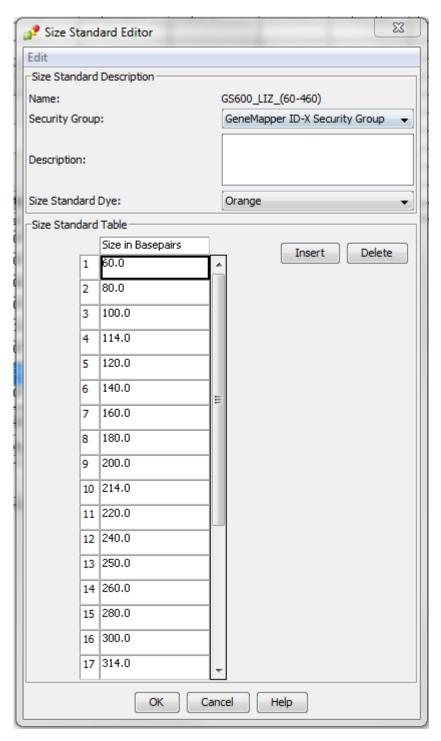
Create a size standard definition file

- 1. Select **Tools GeneMapper ID-X Manager** to open the **GeneMapper ID-X Manager**.
- 2. Click the Size Standards tab, then click New.



- **3.** Specify settings in the **Size Standard Editor**:
 - a. Enter a name as shown in the following figure or enter a new name.
 - **b.** In the **Security Group** field, select the **Security Group** appropriate for your software configuration.
 - c. In the Size Standard Dye field, select Orange.

d. In the **Size Standard Table**, enter the peak sizes that correspond to your size standard.



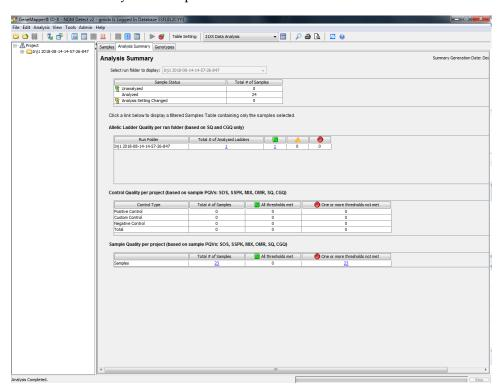
Analyze and edit sample files with GeneMapper ID-X Software

- In the Project window, select Edit > Add Samples to Project, then navigate to the disk or directory that contains the sample files.
- 2. Apply analysis settings to the samples in the project.

Parameter	Settings
Sample Type	Select the sample type.
Analysis Method	Select NGM Detect Analysis Method (or the name of the analysis method you created).
Panel	Select NGM_Detect_Panel.
Size Standard	Use a size range of 60–460 bp for Local Southern size calling method ^[1]

^[1] The NGM Detect[™] kit was originally validated with the GeneScan[™]-600 LIZ[™] Size Standard v2.0. If you use a different size standard, perform the appropriate internal validation studies to support the use of this size standard with the NGM Detect[™] kit.

- **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.
 - The table displays the row of the sample currently being analyzed in green (or red if analysis failed for the sample).
 - The **Analysis Summary** tab is displayed, and the **Genotypes** tab is available when the analysis is complete.



Examine or edit a project

Display electropherogram plots from the Samples and Genotypes tabs of the Project window to examine the data.

For more information on using the GeneMapper $^{\text{\tiny TM}}$ /D-X Software

See "Related documentation" on page 136 for a list of available documents.



Assess the PCR reaction with the Internal Quality Control System

Overview of the Internal Quality Control System	52
Evaluate the PCR reaction	52

Overview of the Internal Quality Control System

The Internal Quality Control System (IQC) is a tool that helps you to assess the PCR reaction and, in conjunction with the STR marker data properties, infer whether a sample is showing signs of degradation or inhibition. The primers for the two IQC markers, IQCS and IQCL, amplify synthetic DNA targets included in the primer mix. IQCS is a low molecular weight amplicon, 70 nt in length. Whereas, IQCL is a high molecular weight amplicon, 456 nt in length.

Note: The allele designation for both IQCS and IQCL will always be 2 (as shown in Figure 2) for all samples. Two alleles for each of the IQC markers are present in the NGM Detect^{TM} Allelic Ladder (designated 1 and 2, see Figure 1). This is because the software dictates that in order for a ladder to be valid, it must have more than one allele per locus.

The IQC system enables you to:

- Confirm the success or failure of the PCR reaction, by looking for the presence or absence of the IQCS and IQCL primer peaks on the electropherogram.
- Determine if PCR inhibitors are present in the PCR reaction, or if PCR reaction conditions are not optimal, by comparing the relative peak heights of IQCS and IQCL.

Evaluate the PCR reaction

To evaluate the PCR performance of the samples, review the relative peak heights of the IQCS and IQCL. Under ideal PCR conditions, the peak height of IQCL is approximately the same or slightly higher than IQCS (termed "balanced" in Table 3). Under sub optimal PCR conditions (for example inhibition), the height of the IQCL is substantially reduced relative to the ICQS. Note that when high inputs of DNA are amplified (greater than 2 ng) some suppression of the IQC peaks may also be seen.

See the following table for outcome scenarios.

Table 3 IQC peak interpretation

Sample DNA profile	IQC peaks	Interpretation	Example	
Balanced	alanced Balanced		See Figure 6	
Ski-slope IQCL peak height significantly decreased or not present		Inhibition	See Figure 7 and Figure 8	
Ski-slope	Balanced	Degraded sample DNA	See Figure 9	
No peaks	Balanced	No DNA or very little sample DNA	See Figure 10	
No peaks	No peaks	PCR failure	_	

Balanced profile

In this example of a balanced profile, the IQC peaks and the DNA profile peaks are balanced. This indicates that PCR has occurred optimally.

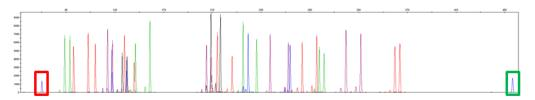


Figure 6 Combined dyes electropherogram showing IQCS and IQCL peaks with 0.5 ng DNA (scaled to 9,000 RFU). The IQCS peak is highlighted by the red box and the IQCL peak is highlighted by the green box.



Ski slope profile with decreased IQCL peak height Figure 7 shows a significantly lower IQCL peak height relative to the IQCS peak height. This indicates that the PCR reaction has been compromised by inhibition. If high levels of PCR inhibition occur, lower IQCL peaks exhibiting some shouldering due to incomplete +A nucleotide addition may be observed. Figure 8 shows the complete absence of an IQCL peak, indicating a high level of inhibition.

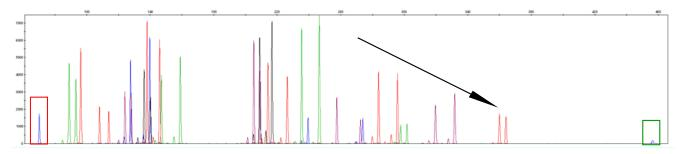


Figure 7 Combined dyes electropherogram for the NGM Detect^{$^{\text{M}}$} kit in the presence of 140 ng/ μ L humic acid. The IQCS peak is highlighted by the red box and the IQCL peak is highlighted by the green box. The arrow indicates the ski slope peak pattern observed in the DNA profile.

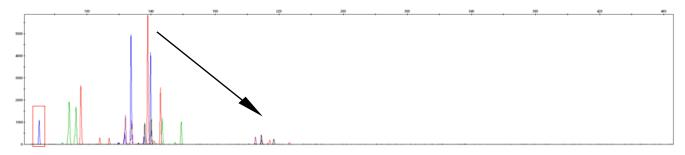


Figure 8 Combined dyes electropherogram for the NGM Detect^{$^{\text{M}}$} kit in the presence of 250 ng/ μ L humic acid. The IQCS peak is highlighted by the red box and the IQCL peak is absent. The arrow indicates the ski slope peak pattern observed in the DNA profile.

Ski slope profile with balanced IQC peaks

The presence of both the IQCS and IQCL with balanced relative peak heights indicates that PCR has occurred optimally.

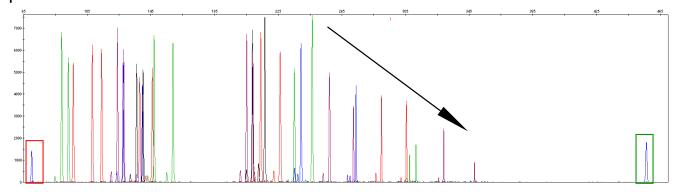


Figure 9 Combined dyes electropherogram from degraded DNA. The IQCS peak is highlighted by the red box and the IQCL peak is highlighted by the green box. The ski slope nature of the DNA profile is highlighted by the arrow.

No sample peaks with balanced IQC peaks

Although there are no DNA profile peaks in the following figure, the presence of both the IQCS and IQCL peaks with balanced relative peak heights indicates that PCR has occurred optimally.

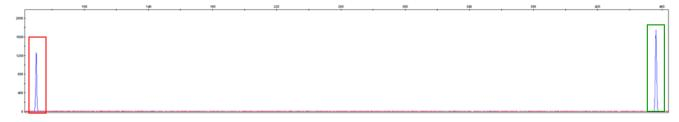


Figure 10 Combined dyes electropherogram showing IQCS and IQCL peaks with 0 ng DNA (scaled to 2,000 RFU). The IQCS peak is highlighted by the red box and the IQCL peak is highlighted by the green box.



Experiments and results

Importance of validation	56
Experiment conditions	56
Laboratory requirements for internal validation	57
Developmental validation	57
Accuracy, precision, and reproducibility	63
Extra peaks in the electropherogram	77
Characterization of loci	88
Species specificity	89
Sensitivity	92
Stability	93
Mixture studies	97
Population data	101

Importance of validation

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

Experiment conditions

We conducted developmental validation experiments according to the updated and revised guidelines from the Scientific Working Group on DNA Analysis Methods (SWGDAM, December 2016). Based on these guidelines, we conducted experiments that comply with guidelines 2.0 and 3.0 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).

We used 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.

Laboratory requirements for internal validation

Each laboratory using this kit must perform internal validation studies. Performance of this kit is supported when used according to the following developmentally validated parameters. Modifications to the protocol should be accompanied by appropriate validation studies performed by the laboratory.

Developmental validation

Except where noted, all developmental validation studies were performed using the GeneAmp[™] PCR System 9700 96-Well thermal cycler or the ProFlex[™] PCR System according to the protocol described in the Perform PCR chapter.

Unless otherwise indicated, the data in this chapter are from the re-validation of the NGM Detect[™] kit after redesigning the IQCL and TH01 markers. The performance of the updated kit formulation assay is fully comparable to that of the original kit. See the *Technical Note: Updated NGM Detect* PCR Amplification Kit: Validation and Comparative Study for studies directly related to soil specificity and a direct comparison of the NGM Detect kit original formulation to the updated formulation.

SWGDAM guideline 2.2.1

"Developmental validation is the acquisition of test data and determination of conditions and limitations of a new or novel DNA methodology for use on forensic, database, known or casework reference samples." (SWGDAM, December 2016)

SWGDAM guideline 3.9.2

"The reaction conditions needed to provide the required degree of specificity and robustness should be determined. These include, but are not limited to, thermal cycling parameters, the concentration of primers, magnesium chloride, DNA polymerase, and other critical reagents." (SWGDAM, December 2016)

SWGDAM guideline 3.9.6

"Criteria for detection of amplified product should be determined based on the platform and/or method." (SWGDAM, December 2016)

PCR components

We examined the concentration of each component of the kit. The concentration of each component was in the range where data indicated that the amplification met the required performance criteria for specificity, sensitivity, and reproducibility. For example, 0.5 ng of DNA Control 007 was amplified in the presence of varying concentrations of magnesium chloride, and the results were analyzed on an Applied Biosystems[™] 3500xL Genetic Analyzer (Figure 11). The performance of the multiplex is most robust within ±20% of the optimal magnesium chloride concentration.

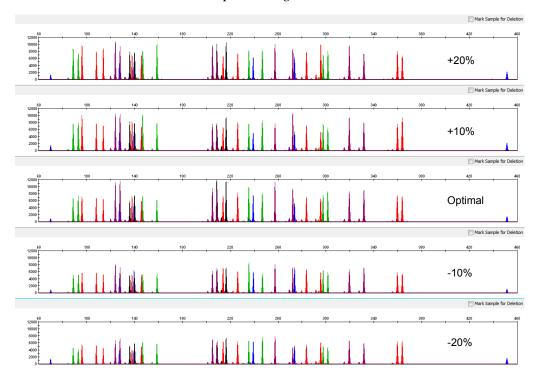


Figure 11 DNA Control 007 (0.5 ng) amplified with the NGM Detect[™] kit in the presence of varying concentrations of magnesium chloride and analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0 to 12,000 RFU).

Thermal cycling temperatures

Thermal cycling parameters were optimized using a Design of Experiments (DOE) approach that seeks to identify the combination of temperatures and hold times that produce the best assay performance. Optimal assay performance was determined through evaluation of several factors, including; evaluation of assay sensitivity, peakheight balance, and resistance to PCR inhibitors.

For example, annealing/extension temperatures of 57, 58, 59, 60, and 61°C were tested using a GeneAmp $^{\text{\tiny M}}$ PCR System 9700 (Figure 12). The PCR products were analyzed using a 3500xL Genetic Analyzer.

6

Of the tested annealing temperatures, 57°C to 61°C produced robust profiles. 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 recommended.

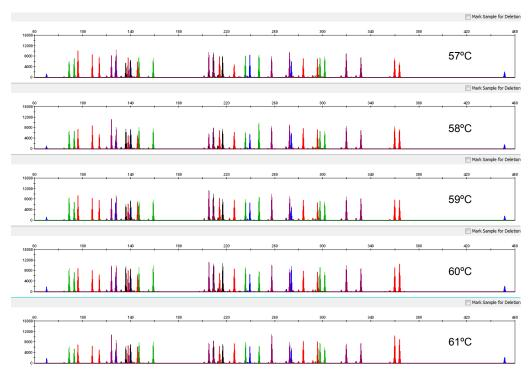


Figure 12 Electropherograms obtained from amplification of 0.5 ng of DNA Control 007 at annealing temperatures of 57, 58, 59, 60, and 61°C, analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0 to 16,000 RFU).

PCR cycle number

Reactions were amplified for 28, 29, 30, 31, and 32 cycles on the GeneAmp[™] PCR System 9700 using 0.5 ng of DNA Control 007. As expected, the amount of PCR product increased with the number of cycles. A full profile was generated for all numbers of thermal cycles (28–32) and off-scale data were collected for several allele peaks at 32 cycles (Figure 13).

We recommend using 30 cycles to optimize signal peak height and minimize artifact or other undesirable peaks.

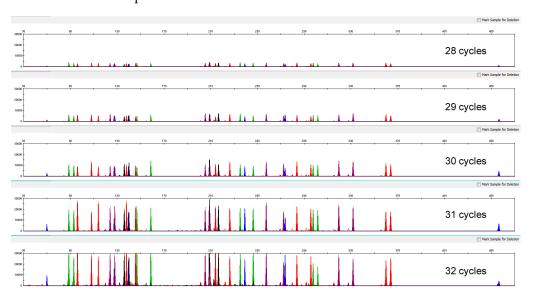


Figure 13 Representative NGM Detect[™] kit profiles obtained from amplification of 0.5 ng of DNA Control 007 using 28, 29, 30, 31, and 32 cycles, analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0 to 30,000 RFU).

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

CE injection time

The effect of injection times on heterozygous peak heights observed for NGM Detect^{$^{\text{TM}}$} assays was studied on three capillary electrophoresis instruments; the 3500 (8-capillary), 3500xL (24-capillary), and 3130xl (16-capillary). As shown in Figure 14 through Figure 16, increasing or decreasing injection times affected profile peak heights in an approximately linear manner. All genomic DNA samples were amplified for 30 cycles.

Our developmental validation studies indicate that the injection conditions that are recorded in Chapter 3 generate profiles from 0.5 ng of input DNA with heterozygous peak height averages between 4,000-10,000 RFU (3500 or 3500xL) or 2,000-4,000 RFU (3130 or 3130xl). This is with no instances of allelic dropout and minimal occurrence of off-scale allele peaks. However, individual CE instrument signal intensities can vary, therefore laboratories are encouraged to optimize injection times. Optimized injection times provide the most efficacious level of assay sensitivity, minimize the

occurrence of off-scale peaks or undesirable artifacts, and do not adversely impact the performance of the size standard and allelic ladder.

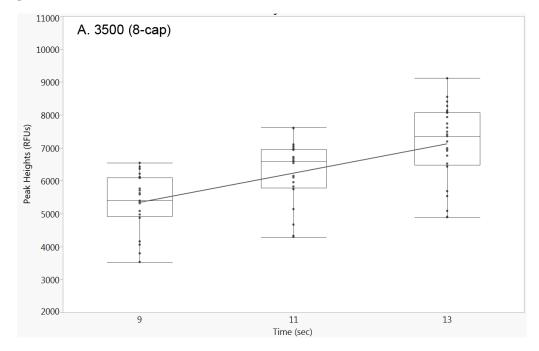


Figure 14 Box plots show the effect of varying injection times (X-axis) on peak heterozygous heights (Y-axis) observed for NGM Detect $^{\text{m}}$ assays with 0.5 ng of human male 007 genomic DNA input on a 3500 Genetic Analyzer.

Note: The data in this figure are from the original formulation of the NGM Detect $^{\text{\tiny TM}}$ kit.

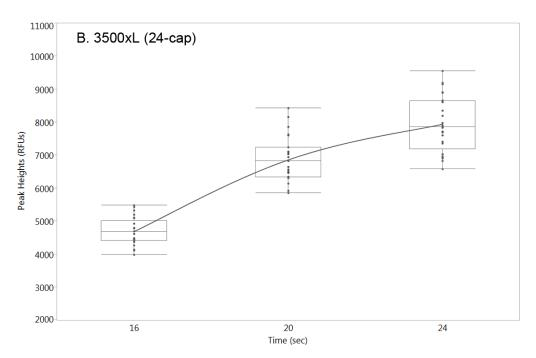


Figure 15 Box plots show the effect of varying injection times (X-axis) on heterozygous peak heights (Y-axis) observed for NGM Detect $^{\text{\tiny M}}$ assays with 0.5 ng of human male 007 genomic DNA input on a 3500xL Genetic Analyzer.

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

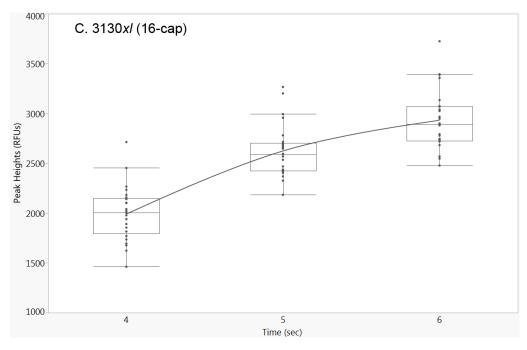


Figure 16 Box plots show the effect of varying injection times (X-axis) on heterozygous peak heights (Y-axis) observed for NGM Detect^{$^{\text{TM}}$} assays with 0.5 ng of human male 007 genomic DNA input on a 3130xl Genetic Analyzer.

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

Accuracy, precision, and reproducibility

SWGDAM guideline 3.5

"Precision and accuracy of the assay should be demonstrated: Precision characterizes the degree of mutual agreement among a series of individual measurements, values and/or results. Precision depends only on the distribution of random errors and does not relate to the true value or specified value. The measure of precision is usually expressed in terms of imprecision and computed as a standard deviation of the test results. Accuracy is the degree of conformity of a measured quantity to its actual (true) value. Accuracy of a measuring instrument is the ability of a measuring instrument to give responses close to a true value." (SWGDAM, December 2016)

Accuracy observation

Laser-induced fluorescence detection of length polymorphism at short tandem repeat loci is not a novel methodology (Holt *et al.*, 2000; Wallin *et al.*, 2002). However, accuracy and reproducibility of profiles have been determined from various sample types.

Figure 17, Figure 18, and Figure 19 show the size differences that were observed between sample alleles and allelic ladder alleles on the Applied Biosystems $^{\text{\tiny TM}}$ 3130 xl, 3500, and 3500xL Genetic Analyzers with POP-4 $^{\text{\tiny TM}}$ Polymer. The X-axis in the following figures represents the nominal nucleotide sizes for the NGM Detect $^{\text{\tiny TM}}$ 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, irrespective of the capillary electrophoresis platforms.

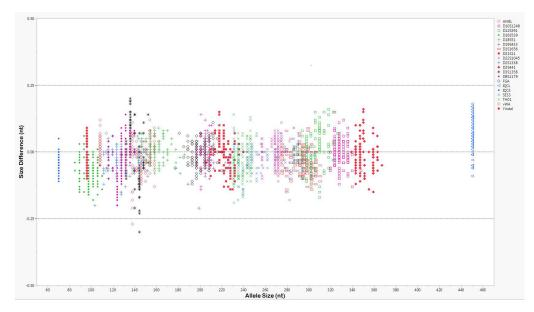


Figure 17 Allele size vs. Allelic Ladder sizing for 81 samples analyzed on a 3130 xl Genetic Analyzer. Size and ladder sizing for the NGM Detect kit were calculated using the GeneScan $^{\text{TM}}$ –600 LIZ Size Standard v2.0.

Note: Except for the TH01 marker data, the data in this figure are from the original formulation of the NGM $Detect^{\mathsf{TM}}$ kit.

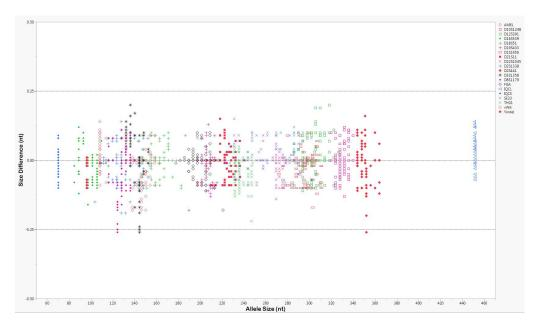


Figure 18 Allele size vs. Allelic Ladder sizing for 81 samples analyzed on a 3500 Genetic Analyzer. Size and ladder sizing for the NGM Detect $^{\text{\tiny TM}}$ kit were calculated using the GeneScan $^{\text{\tiny TM}}$ -600 LIZ $^{\text{\tiny TM}}$ Size Standard v2.0.

Note: Except for the TH01 marker data, the data in this figure are from the original formulation of the NGM $Detect^{\mathsf{TM}}$ kit.

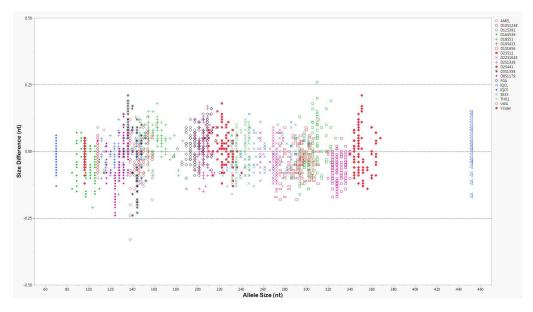


Figure 19 Allele size vs. Allelic Ladder sizing for 81 samples analyzed on a 3500xL Genetic Analyzer. Size and ladder sizing for the NGM Detect $^{\text{\tiny TM}}$ kit were calculated using the GeneScan $^{\text{\tiny TM}}$ –600 LIZ $^{\text{\tiny TM}}$ Size Standard v2.0.

Note: Except for the TH01 marker data, the data in this figure are from the original formulation of the NGM $Detect^{T}$ kit.

Precision and size window description

Sizing precision enables the determination of accurate and reliable genotypes. The recommended method for genotyping is to use a ± 0.5 -nt "window" around the size obtained for each allele in the 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 either:

- An "off-ladder" allele, that is, an allele of a size that is not represented in the allelic ladder.
- An allele that does correspond to an allele in the allelic ladder, 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.

Precision and size window observation

Table 4 lists typical precision results obtained from multiple runs of the NGM Detect Allelic Ladder using the GeneScan^{\mathbb{T}}-600 LIZ^{\mathbb{T}} Size Standard v2.0. The results were obtained within a set of injections on a single capillary array. The number of repeated injections for each genetic analyzer platform is shown in the following table:

CE platform	Capillaries	# Injections	Sizing method
3130 <i>xl</i>	16/injection	5	Local Southern, 60–460 bp
3500	8/injection	12	Local Southern, 60–460 bp
3500xL	24/injection	4	Local Southern, 60–460 bp

The mean sizes and the standard deviation for the allele sizing were calculated for all the alleles in each run (Table 4). The mean range and the standard deviation range show the lowest and highest values obtained across multiple runs.

Sample alleles can 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 with the smallest standard deviations in sizing. The figures in "Accuracy observation" on page 63 illustrate the tight clustering of allele sizes obtained on the Applied Biosystems genetic analyzers, 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 to an allele in the Allelic Ladder. Repeat analysis, when necessary, provides an added level of confidence in the final allele assignment.

GeneMapper $^{\text{\tiny TM}}$ *ID-X* Software automatically flags sample alleles that do not size within the prescribed window around an allelic ladder allele by labeling 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 several 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.

IMPORTANT! To minimize the variation in sizing between runs and to ensure accurate genotyping, follow the guidelines in "Allelic ladder requirements for data analysis" on page 34 and use allelic ladders obtained from the same run as samples to analyze the samples.

For more information on precision and genotyping, see (Lazaruk *et al.*, 1998; Mansfield *et al.*, 1998).

Note: The IQCS and IQCL markers were omitted from this study because they are not used for genotyping.

Table 4 Precision results of multiple runs of the NGM Detect[™] Allelic Ladder

Note: Except for the TH01 marker data, the data in this table are from the original formulation of the NGM Detect $^{\text{TM}}$ kit.

	To initiation of the Now Beteet - Kit.						
	313	0 <i>xl</i>	35	00	350	0xL	
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation	
D2S133	8						
11	90.76-90.87	0.023-0.034	91.05-91.13	0.003-0.040	90.96-91.03	0.028-0.035	
12	94.90-95.00	0.021-0.031	95.17-95.24	0.021-0.041	95.10-95.16	0.024-0.034	
13	99.01-99.13	0.023-0.035	99.31-99.38	0.001-0.041	99.22-99.28	0.027-0.034	
14	103.16-103.28	0.020-0.036	103.45-103.52	0.028-0.044	103.38-103.44	0.031-0.038	
15	107.09-107.19	0.025-0.037	107.31-107.38	0.015-0.058	107.23-107.28	0.028-0.037	
16	111.29-111.38	0.028-0.033	111.55-111.60	0.005-0.043	111.45-111.50	0.025-0.041	
17	115.34-115.43	0.022-0.035	115.56–115.62	0.001-0.032	115.49-115.52	0.011-0.031	
18	119.29-119.36	0.012-0.033	119.52-119.57	0.024-0.042	119.43-119.47	0.015-0.040	
19	123.27-123.36	0.020-0.034	123.50-123.56	0.004-0.041	123.42-123.47	0.034-0.036	
20	127.26-127.35	0.026-0.041	127.51-127.57	0.019-0.054	127.41-127.46	0.028-0.042	
21	131.28-131.39	0.024-0.039	131.53-131.62	0.017-0.051	131.46-131.51	0.029-0.039	
22	135.34-135.44	0.023-0.038	135.56-135.63	0.011-0.046	135.50-135.53	0.027-0.035	
23	139.39-139.49	0.020-0.040	139.62-139.69	0.001-0.048	139.54-139.59	0.002-0.033	
24	143.51-143.61	0.020-0.040	143.77-143.82	0.008-0.043	143.68-143.74	0.026-0.042	
25	147.63-147.74	0.017-0.031	147.89-147.94	0.016-0.048	147.79–147.85	0.024-0.043	
26	151.76-151.85	0.023-0.029	152.02-152.06	0.018-0.041	151.92-151.97	0.030-0.036	
27	155.91–156.01	0.018-0.031	156.15-156.20	0.028-0.043	156.07-156.11	0.026-0.040	

	313	80 <i>xl</i>	35	00	350	0xL
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
28	160.45-160.49	0.008-0.033	160.51-160.57	0.007-0.042	160.45-160.47	0.037-0.040
SE33						
4.2	188.38-188.44	0.024-0.037	188.22-188.26	0.018-0.044	188.17–188.18	0.034-0.036
6.3	197.70-197.76	0.022-0.038	197.54–197.6	0.005-0.045	197.49-197.50	0.027-0.038
8	202.76-202.80	0.022-0.032	202.58-202.63	0.011-0.043	202.53-202.55	0.022-0.033
9	206.78-206.85	0.023-0.040	206.60-206.65	0.017-0.044	206.56-206.56	0.030-0.036
11	214.87-214.92	0.031-0.036	214.65-214.69	0.035-0.051	214.59-214.62	0.029-0.038
12	219.05-219.12	0.021-0.037	218.83-218.88	0.033-0.053	218.78-218.80	0.035-0.042
13	223.16-223.20	0.032-0.047	222.92-222.97	0.021-0.050	222.88-222.90	0.033-0.043
14	227.22-227.29	0.032-0.046	227.01-227.06	0.011-0.054	226.96-226.98	0.033-0.049
15	231.26-231.34	0.028-0.047	231.06-231.09	0.018-0.070	231.01-231.03	0.037-0.045
16	235.40-235.46	0.024-0.041	235.16-235.22	0.013-0.053	235.14-235.14	0.039-0.044
17	239.50-239.56	0.029-0.045	239.27-239.32	0.031-0.048	239.23-239.24	0.039-0.045
18	243.65-243.71	0.020-0.031	243.38-243.43	0.004-0.044	243.35-243.36	0.027-0.042
19	247.78-247.84	0.030-0.037	247.47-247.54	0.009-0.049	247.47-247.48	0.026-0.035
20	251.90-251.95	0.026-0.036	251.58-251.65	0.009-0.049	251.57-251.59	0.033-0.037
20.2	253.86-253.92	0.021-0.041	253.56-253.63	0.025-0.045	253.55-253.56	0.028-0.036
21	255.86-255.93	0.028-0.038	255.57-255.63	0.015-0.048	255.55-255.57	0.028-0.040
21.2	257.83-257.90	0.026-0.042	257.55-257.60	0.031-0.043	257.52-257.54	0.021-0.030
22.2	261.81-261.88	0.019-0.043	261.53-261.58	0.005-0.052	261.51-261.53	0.030-0.039
23.2	265.91-265.98	0.025-0.047	265.62-265.68	0.027-0.055	265.60-265.62	0.030-0.036
24.2	269.97-270.02	0.026-0.041	269.66-269.72	0.025-0.055	269.64-269.66	0.035-0.041
25.2	273.95-274.00	0.025-0.036	273.65-273.69	0.034-0.053	273.62-273.64	0.031-0.045
26.2	278.10-278.16	0.029-0.043	277.8-277.84	0.007-0.058	277.78-277.80	0.033-0.038
27.2	282.16-282.22	0.027-0.047	281.84-281.89	0.005-0.058	281.81-281.84	0.026-0.050
28.2	286.14-286.21	0.022-0.045	285.83-285.86	0.023-0.054	285.80-285.82	0.027-0.038
29.2	290.15-290.22	0.028-0.046	289.85-289.89	0.016-0.052	289.82-289.83	0.031-0.036
30.2	294.20-294.24	0.028-0.039	293.87-293.92	0.029-0.066	293.85-293.85	0.037-0.045
31.2	298.16-298.24	0.028-0.045	297.84-297.92	0.001-0.050	297.82-297.83	0.032-0.043

	313	0 <i>xl</i>	35	00	350	0xL
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
32.2	302.17-302.23	0.027-0.041	301.83-301.90	0.004-0.051	301.81-301.83	0.042-0.045
33.2	306.14-306.20	0.025-0.050	305.81-305.86	0.007-0.043	305.76-305.80	0.030-0.033
34.2	310.16-310.22	0.033-0.045	309.81-309.88	0.007-0.048	309.78-309.80	0.034-0.043
35	312.20-312.27	0.025-0.050	311.84-311.89	0.005-0.047	311.82-311.83	0.034-0.036
35.2	314.24-314.30	0.027-0.041	313.87-313.94	0.001-0.050	313.83-313.87	0.037-0.048
36	316.37-316.43	0.024-0.042	316-316.04	0.010-0.038	315.97-315.99	0.024-0.048
37	320.60-320.66	0.032-0.048	320.21-320.27	0.035-0.053	320.19-320.23	0.037-0.048
38	324.76-324.83	0.035-0.054	324.35-324.41	0.009-0.064	324.35-324.38	0.042-0.048
39	328.85-328.92	0.027-0.047	328.45-328.51	0.029-0.055	328.45-328.48	0.029-0.039
42	341.00-341.06	0.031-0.048	340.58-340.65	0.001-0.064	340.57-340.59	0.037-0.044
D16S53	19					
5	72.32-72.38	0.032-0.041	72.00-72.05	0.009-0.051	71.93-71.98	0.024-0.031
8	84.95-84.99	0.024-0.040	84.64-84.70	0.008-0.044	84.58-84.61	0.027-0.040
9	89.15-89.18	0.020-0.034	88.84-88.9	0.004-0.045	88.78-88.82	0.027-0.041
10	93.30-93.33	0.022-0.036	92.99-93.05	0.004-0.040	92.94-92.97	0.028-0.040
11	97.45-97.49	0.020-0.039	97.15-97.22	0.007-0.045	97.11-97.14	0.025-0.035
12	101.63-101.67	0.026-0.033	101.33-101.39	0.030-0.046	101.28-101.33	0.034-0.039
13	105.83-105.86	0.021-0.043	105.53-105.59	0.009-0.047	105.47-105.52	0.038-0.039
14	109.97-110.01	0.027-0.042	109.66-109.74	0.032-0.048	109.63-109.67	0.030-0.043
15	114.08-114.11	0.024-0.036	113.77-113.85	0.028-0.051	113.72-113.78	0.027-0.038
D18S51						
7	127.11-127.18	0.030-0.038	127.52-127.58	0.014-0.046	127.44-127.47	0.028-0.042
9	134.69-134.79	0.022-0.047	135.21-135.27	0.010-0.051	135.12-135.17	0.028-0.045
10	138.52-138.63	0.025-0.051	139.07-139.15	0.001-0.041	139.00-139.06	0.037-0.042
10.2	140.66-140.77	0.028-0.054	141.21-141.30	0.001-0.042	141.15-141.21	0.035-0.043
11	142.41-142.53	0.023-0.047	143.00-143.10	0.008-0.052	142.93-142.99	0.020-0.038
12	146.33-146.45	0.023-0.050	146.95–147.03	0.020-0.044	146.88-146.95	0.029-0.038
13	150.27-150.39	0.027-0.042	150.90-151.01	0.024-0.045	150.83-150.91	0.031-0.038
13.2	152.41-152.53	0.027-0.050	153.05-153.16	0.013-0.046	153.00-153.07	0.033-0.044

	313	i0xl	35	00	3500xL	
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
14	154.21–154.33	0.030-0.048	154.86-154.96	0.024-0.046	154.80-154.87	0.032-0.036
14.2	156.35-156.48	0.021-0.040	157.02-157.14	0.004-0.047	156.96-157.03	0.035-0.043
15	158.15-158.27	0.032-0.051	158.83-158.94	0.001-0.043	158.76-158.83	0.040-0.045
16	162.05-162.18	0.027-0.049	162.76-162.86	0.004-0.043	162.69-162.77	0.030-0.037
17	165.92-166.07	0.027-0.054	166.64-166.77	0.003-0.039	166.58-166.66	0.029-0.046
18	169.80-169.93	0.034-0.049	170.54-170.66	0.016-0.044	170.49-170.57	0.030-0.039
19	173.66-173.81	0.031-0.049	174.44-174.55	0.004-0.044	174.39-174.47	0.029-0.045
20	177.54-177.69	0.035-0.051	178.34–178.48	0.003-0.047	178.28-178.36	0.032-0.040
21	181.47-181.63	0.039-0.049	182.30-182.45	0.006-0.044	182.26-182.35	0.035-0.044
22	185.33–185.52	0.035-0.047	186.21–186.37	0.013-0.042	186.17–186.26	0.033-0.046
23	189.25-189.44	0.030-0.067	190.17-190.32	0.022-0.046	190.12-190.22	0.032-0.042
24	193.19-193.37	0.032-0.054	194.13-194.26	0.021-0.050	194.07–194.17	0.029-0.055
25	197.14-197.32	0.030-0.064	198.11-198.25	0.005-0.048	198.05-198.15	0.033-0.047
26	200.98-201.15	0.038-0.065	201.92-202.07	0.024-0.048	201.88-201.98	0.034-0.043
27	204.81-204.98	0.035-0.067	205.76-205.91	0.020-0.040	205.72-205.83	0.038-0.046
TH01						
4	222.50-222.65	0.035-0.073	223.26-223.39	0.004-0.049	223.19-223.31	0.035-0.041
5	226.48-226.64	0.029-0.076	227.24-227.39	0.011-0.060	227.19-227.32	0.031-0.040
6	230.47-230.65	0.034-0.057	231.25-231.39	0.029-0.049	231.19-231.31	0.027-0.044
7	234.46-234.65	0.028-0.059	235.27-235.39	0.008-0.048	235.18-235.31	0.033-0.041
8	238.47-238.65	0.033-0.068	239.25-239.40	0.005-0.050	239.19-239.32	0.029-0.043
9	242.55-242.71	0.028-0.065	243.34-243.47	0.027-0.053	243.26-243.38	0.025-0.040
9.3	245.61-245.78	0.035-0.058	246.38-246.51	0.005-0.049	246.31-246.43	0.034-0.047
10	246.63-246.81	0.033-0.074	247.40-247.54	0.024-0.050	247.33-247.45	0.031-0.041
11	250.70-250.86	0.036-0.066	251.43-251.54	0.003-0.046	251.36-251.47	0.034-0.043
12	254.62-254.79	0.031-0.069	255.36-255.48	0.016-0.046	255.29-255.40	0.030-0.044
13	258.53-258.71	0.036-0.063	259.28-259.39	0.003-0.045	259.19-259.31	0.033-0.040
13.3	261.46-261.65	0.042-0.065	262.21-262.34	0.005-0.038	262.15-262.27	0.027-0.042

	3130 <i>xl</i>		3500		3500xL	
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
D12539	1					
14	281.70-281.76	0.022-0.037	281.71-281.75	0.039-0.054	281.68-281.70	0.033-0.041
15	285.72-285.76	0.026-0.033	285.73-285.75	0.010-0.048	285.69-285.71	0.030-0.043
16	289.68-289.72	0.023-0.040	289.68-289.71	0.001-0.049	289.65-289.67	0.030-0.038
17	293.71-293.77	0.018-0.037	293.71-293.76	0.001-0.046	293.68-293.69	0.027-0.041
18	297.70-297.75	0.026-0.040	297.7-297.74	0.001-0.050	297.67-297.69	0.036-0.040
19	301.69-301.74	0.020-0.042	301.70-301.75	0.024-0.048	301.65-301.67	0.032-0.040
19.3	304.64-304.68	0.021-0.033	304.62-304.67	0.011-0.047	304.58-304.61	0.031-0.037
20	305.62-305.66	0.025-0.037	305.61-305.65	0.007-0.048	305.56-305.59	0.030-0.039
21	309.65-309.70	0.022-0.043	309.65-309.71	0.012-0.054	309.60-309.64	0.036-0.044
22	313.76-313.81	0.028-0.041	313.77-313.79	0.031-0.050	313.71-313.74	0.034-0.040
23	318.00-318.04	0.022-0.037	317.97-318.00	0.032-0.051	317.92-317.96	0.026-0.039
24	322.19-322.22	0.026-0.046	322.13-322.18	0.004-0.056	322.11-322.13	0.016-0.047
25	326.3-326.33	0.022-0.041	326.24-326.28	0.011-0.070	326.21-326.23	0.034-0.052
26	330.39-330.42	0.027-0.033	330.34-330.38	0.029-0.067	330.31-330.34	0.037-0.048
27	334.42-334.46	0.025-0.039	334.38-334.40	0.016-0.060	334.34-334.38	0.027-0.042
D3S135	8					
9	112.46-112.49	0.019-0.047	112.54-112.59	0.038-0.046	112.50-112.53	0.017-0.035
10	116.39-116.42	0.021-0.027	116.47-116.51	0.001-0.036	116.43-116.45	0.024-0.032
11	120.26-120.29	0.018-0.035	120.33-120.37	0.028-0.044	120.29-120.30	0.022-0.037
12	124.05-124.08	0.023-0.034	124.11-124.15	0.007-0.049	124.08-124.09	0.023-0.044
13	128.11-128.15	0.024-0.038	128.17-128.22	0.025-0.049	128.14-128.15	0.031-0.037
14	132.04-132.09	0.016-0.035	132.12-132.16	0.026-0.050	132.09-132.10	0.029-0.038
15	135.93-135.96	0.024-0.034	136.01-136.04	0.006-0.037	135.98-135.98	0.030-0.035
16	140.07-140.11	0.017-0.035	140.16-140.20	0.001-0.046	140.13-140.13	0.039-0.043
17	144.27–144.29	0.020-0.036	144.33-144.37	0.034-0.052	144.31-144.32	0.027-0.037
18	148.33–148.36	0.017-0.031	148.4-148.44	0.029-0.049	148.38-148.38	0.028-0.033
19	152.35–152.37	0.019-0.030	152.40-152.45	0.027-0.045	152.38-152.39	0.032-0.039
20	156.66–156.70	0.023-0.037	156.70–156.76	0.008-0.038	156.69-156.71	0.020-0.037

Allele	3130 <i>xl</i>		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
FGA						
13	165.50-165.58	0.029-0.039	165.95–166.04	0.015-0.047	165.93–165.97	0.023-0.040
14	169.35-169.45	0.030-0.036	169.84-169.91	0.003-0.043	169.81-169.86	0.030-0.035
15	173.21-173.31	0.031-0.039	173.72-173.83	0.009-0.043	173.70-173.74	0.028-0.038
16	177.06-177.17	0.032-0.041	177.59-177.71	0.004-0.047	177.58-177.63	0.031-0.041
17	180.94-181.05	0.026-0.042	181.51-181.61	0.005-0.047	181.48-181.55	0.027-0.047
18	184.85–184.97	0.032-0.045	185.45-185.56	0.013-0.051	185.43-185.50	0.030-0.039
19	188.76–188.89	0.024-0.040	189.38-189.51	0.021-0.039	189.38-189.45	0.031-0.042
20	192.67-192.80	0.032-0.042	193.34-193.44	0.010-0.049	193.31-193.39	0.028-0.049
21	196.56–196.70	0.030-0.050	197.23-197.38	0.004-0.046	197.24-197.32	0.035-0.043
22	200.46-200.59	0.034-0.054	201.15-201.27	0.027-0.054	201.13-201.22	0.026-0.043
23	204.28-204.39	0.033-0.049	204.97-205.11	0.020-0.044	204.96-205.05	0.031-0.050
24	208.09-208.22	0.037-0.048	208.83-208.93	0.011-0.056	208.81-208.89	0.035-0.048
25	211.94-212.05	0.027-0.048	212.67-212.82	0.004-0.048	212.67-212.74	0.039-0.050
26	215.83-215.95	0.029-0.061	216.58-216.74	0.023-0.041	216.59-216.69	0.037-0.043
27	219.83-219.97	0.034-0.059	220.62-220.79	0.001-0.050	220.64-220.73	0.036-0.050
28	223.71-223.86	0.035-0.058	224.55-224.68	0.011-0.053	224.54-224.64	0.037-0.040
29	227.55-227.70	0.034-0.063	228.41-228.58	0.016-0.061	228.43-228.53	0.030-0.056
30	231.34-231.50	0.034-0.056	232.23-232.40	0.015-0.057	232.25-232.36	0.035-0.056
30.2	233.71-233.86	0.030-0.054	234.60-234.76	0.014-0.050	234.62-234.73	0.039-0.052
31.2	237.59-237.74	0.029-0.062	238.50-238.68	0.003-0.050	238.52-238.64	0.040-0.051
32.2	241.50-241.66	0.038-0.053	242.46-242.62	0.034-0.050	242.47-242.59	0.038-0.047
33.2	245.48-245.64	0.035-0.049	246.43-246.61	0.022-0.053	246.47-246.56	0.041-0.052
42.2	280.49-280.68	0.034-0.067	281.59-281.80	0.001-0.048	281.61-281.75	0.038-0.052
43.2	284.32-284.52	0.041-0.073	285.46-285.67	0.016-0.047	285.47-285.62	0.032-0.052
44.2	288.33-288.51	0.041-0.070	289.46-289.66	0.016-0.053	289.48-289.62	0.033-0.055
45.2	292.23-292.41	0.040-0.070	293.35-293.57	0.001-0.055	293.38-293.52	0.033-0.048
46.2	295.73-295.93	0.037-0.065	296.95-297.17	0.004-0.051	296.98-297.12	0.038-0.059
47.2	299.50-299.7	0.037-0.073	300.72-300.95	0.007-0.059	300.76-300.90	0.036-0.065

Allele	3130 <i>xl</i>		3500		3500xL				
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation			
48.2	303.40-303.60	0.034-0.075	304.64-304.87	0.018-0.053	304.67-304.82	0.040-0.055			
50.2	311.14-311.37	0.048-0.081	312.47-312.69	0.007-0.063	312.46-312.64	0.043-0.057			
51.2	314.87-315.12	0.052-0.083	316.30-316.56	0.021-0.066	316.34-316.51	0.039-0.055			
Yindel									
1	90.97-91.00	0.021-0.031	91.04-91.10	0.028-0.044	90.97-90.98	0.030-0.033			
2	96.09-96.12	0.022-0.033	96.18-96.22	0.001-0.035	96.10-96.11	0.022-0.035			
AMEL									
Х	108.18-108.22	0.024-0.039	108.08-108.16	0.007-0.046	108.05-108.09	0.030-0.035			
Υ	113.96-114.01	0.001-0.035	113.94-113.99	0.028-0.042	113.91-113.94	0.025-0.035			
vWA									
11	125.73-125.77	0.017-0.040	125.79-125.85	0.015-0.046	125.74-125.78	0.024-0.038			
12	129.63-129.66	0.021-0.036	129.72-129.76	0.026-0.046	129.67-129.69	0.026-0.036			
13	133.62-133.65	0.023-0.039	133.74-133.75	0.017-0.050	133.67-133.70	0.031-0.036			
14	137.82-137.86	0.021-0.035	137.96-137.99	0.004-0.045	137.92-137.93	0.025-0.040			
15	141.72-141.75	0.024-0.036	141.87-141.90	0.005-0.045	141.81-141.83	0.025-0.030			
16	145.81-145.83	0.022-0.035	145.96-146.01	0.026-0.043	145.92-145.93	0.027-0.036			
17	149.88-149.92	0.024-0.035	150.05-150.10	0.028-0.042	150.02-150.03	0.034-0.040			
18	153.94-153.97	0.027-0.035	154.12-154.16	0.021-0.046	154.07-154.08	0.028-0.040			
19	158.03-158.07	0.023-0.040	158.23-158.27	0.004-0.046	158.18-158.20	0.031-0.038			
20	162.08-162.13	0.022-0.048	162.29-162.32	0.024-0.045	162.24-162.27	0.023-0.037			
21	166.07-166.11	0.024-0.033	166.26-166.33	0.026-0.045	166.24-166.25	0.022-0.038			
22	170.08-170.11	0.016-0.036	170.29-170.34	0.020-0.041	170.25-170.27	0.032-0.035			
23	173.97-174.03	0.020-0.032	174.20-174.26	0.010-0.047	174.17-174.18	0.031-0.039			
24	178.37-178.42	0.029-0.037	178.59-178.64	0.003-0.046	178.56-178.57	0.029-0.038			
D21S11									
24	197.75-197.85	0.024-0.039	198.20-198.26	0.032-0.046	198.16-198.20	0.034-0.045			
24.2	199.77-199.87	0.033-0.040	200.23-200.30	0.001-0.046	200.19-200.24	0.023-0.042			
25	201.76-201.85	0.023-0.045	202.20-202.25	0.031-0.043	202.16-202.2	0.027-0.035			
26	205.74-205.81	0.020-0.046	206.16-206.22	0.024-0.050	206.12-206.16	0.031-0.044			

	313	0xl	3500		3500xL	
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
27	209.74-209.81	0.025-0.043	210.16-210.23	0.031-0.047	210.12-210.17	0.037-0.046
28	213.69-213.75	0.025-0.037	214.11-214.16	0.031-0.048	214.08-214.11	0.025-0.042
28.2	215.69-215.75	0.029-0.044	216.11-216.19	0.025-0.046	216.09-216.13	0.028-0.040
29	217.69-217.77	0.024-0.032	218.13-218.22	0.025-0.045	218.13-218.16	0.029-0.042
29.2	219.81-219.88	0.027-0.045	220.23-220.31	0.001-0.049	220.22-220.26	0.030-0.041
30	221.77-221.86	0.024-0.037	222.23-222.3	0.005-0.043	222.21-222.25	0.032-0.047
30.2	223.77-223.84	0.027-0.045	224.20-224.29	0.008-0.056	224.20-224.24	0.034-0.045
31	225.78-225.86	0.027-0.049	226.26-226.32	0.021-0.059	226.24-226.28	0.038-0.043
31.2	227.76-227.85	0.030-0.041	228.23-228.32	0.011-0.054	228.21-228.28	0.037-0.039
32	229.78-229.86	0.033-0.047	230.28-230.35	0.022-0.056	230.26-230.31	0.038-0.043
32.2	231.76-231.86	0.030-0.049	232.27-232.34	0.019-0.055	232.25-232.29	0.027-0.042
33	233.79-233.89	0.027-0.043	234.28-234.38	0.009-0.048	234.28-234.33	0.036-0.038
33.2	235.74-235.83	0.031-0.043	236.27-236.33	0.011-0.046	236.23-236.29	0.037-0.040
34	237.87-237.98	0.023-0.040	238.38-238.47	0.007-0.048	238.38-238.42	0.017-0.043
34.2	239.80-239.89	0.036-0.041	240.33-240.40	0.001-0.048	240.30-240.37	0.033-0.044
35	241.92-242.02	0.029-0.044	242.46-242.54	0.030-0.052	242.43-242.49	0.036-0.041
35.2	243.92-244.00	0.029-0.038	244.43-244.51	0.024-0.053	244.41-244.47	0.028-0.041
36	245.94-246.05	0.025-0.037	246.46-246.55	0.004-0.048	246.46-246.49	0.033-0.040
37	250.09-250.18	0.032-0.039	250.59-250.67	0.001-0.046	250.58-250.63	0.028-0.040
38	253.95-254.05	0.024-0.044	254.46-254.53	0.018-0.048	254.45-254.51	0.028-0.043
D1S165	6					
9	267.24-267.33	0.029-0.037	267.69-267.75	0.019-0.049	267.64-267.68	0.035-0.040
10	271.29-271.37	0.027-0.034	271.71-271.79	0.021-0.047	271.67-271.72	0.035-0.037
11	275.31-275.40	0.026-0.041	275.76-275.82	0.010-0.048	275.70-275.76	0.031-0.040
12	279.35-279.42	0.021-0.039	279.76-279.83	0.001-0.049	279.71-279.76	0.035-0.046
13	283.38-283.45	0.017-0.043	283.80-283.86	0.001-0.051	283.73-283.79	0.031-0.039
14	287.35-287.40	0.022-0.035	287.75-287.82	0.005-0.037	287.70-287.74	0.030-0.041
14.3	290.42-290.48	0.027-0.040	290.80-290.86	0.013-0.055	290.75-290.81	0.038-0.047
15	291.31-291.37	0.018-0.037	291.72-291.79	0.005-0.050	291.67-291.72	0.031-0.040

	313	0 <i>xl</i>	35	00	350	0xL
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
15.3	294.41-294.45	0.025-0.037	294.80-294.85	0.010-0.050	294.74-294.78	0.035-0.047
16	295.31-295.38	0.020-0.038	295.75-295.80	0.008-0.048	295.66-295.72	0.033-0.039
16.3	298.41-298.45	0.029-0.036	298.78-298.84	0.031-0.049	298.71-298.76	0.038-0.045
17	299.24-299.29	0.022-0.043	299.64-299.72	0.001-0.052	299.56-299.61	0.018-0.044
17.3	302.34-302.39	0.028-0.040	302.71-302.77	0.023-0.052	302.64-302.69	0.023-0.042
18	303.19-303.24	0.021-0.037	303.62-303.65	0.007-0.047	303.53-303.58	0.030-0.036
18.3	306.24-306.32	0.024-0.040	306.66-306.70	0.018-0.048	306.57-306.64	0.026-0.048
19.3	310.25-310.32	0.026-0.044	310.65-310.70	0.007-0.057	310.57-310.63	0.029-0.036
20.3	314.22-314.30	0.008-0.041	314.65-314.71	0.001-0.048	314.56-314.62	0.018-0.045
D2S441						
8	334.96-335.09	0.033-0.047	335.78-335.91	0.031-0.050	335.70-335.84	0.027-0.045
9	338.84-338.99	0.035-0.049	339.68-339.78	0.031-0.053	339.57-339.71	0.018-0.053
10	342.87-343.01	0.029-0.049	343.71-343.87	0.032-0.054	343.64-343.78	0.035-0.047
11	346.91-347.05	0.035-0.046	347.80-347.91	0.016-0.056	347.68-347.82	0.030-0.049
11.3	349.92-350.06	0.032-0.061	350.83-350.96	0.016-0.049	350.72-350.87	0.026-0.054
12	351.00-351.15	0.026-0.054	351.91-352.03	0.021-0.055	351.80-351.94	0.035-0.053
13	354.97-355.12	0.032-0.051	355.88-356.01	0.010-0.050	355.78-355.94	0.027-0.047
14	358.95-359.10	0.036-0.056	359.88-360	0.001-0.053	359.78-359.90	0.037-0.048
15	362.94-363.09	0.040-0.058	363.86-363.99	0.010-0.051	363.77-363.92	0.040-0.044
16	366.93-367.07	0.037-0.049	367.83-367.97	0.024-0.051	367.77-367.91	0.033-0.039
17	370.94-371.08	0.030-0.051	371.86-372.01	0.035-0.049	371.79-371.94	0.032-0.043
D8S117	9					
5	95.58-95.63	0.025-0.030	95.79-95.84	0.027-0.045	95.73-95.76	0.025-0.035
6	99.66-99.71	0.016-0.032	99.88-99.92	0.001-0.042	99.83-99.84	0.016-0.034
7	103.81-103.86	0.023-0.033	104.04-104.09	0.010-0.047	103.98-104.01	0.028-0.036
8	107.91-107.96	0.023-0.037	108.15-108.19	0.007-0.044	108.09-108.10	0.033-0.042
9	111.97-112.02	0.017-0.034	112.21-112.26	0.007-0.045	112.17-112.18	0.025-0.035
10	115.96-116.00	0.017-0.033	116.19-116.24	0.001-0.039	116.15-116.16	0.019-0.026
11	119.87-119.91	0.024-0.031	120.12-120.16	0.001-0.051	120.07-120.08	0.014-0.024

	313	80 <i>xl</i>	35	00	350	0xL
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
12	123.91-123.96	0.027-0.033	124.17-124.22	0.007-0.052	124.13-124.15	0.026-0.040
13	127.72-127.79	0.022-0.038	128.01-128.06	0.021-0.043	127.96-127.98	0.030-0.041
14	131.64-131.70	0.021-0.038	131.94-132.00	0.019-0.050	131.91-131.93	0.031-0.041
15	135.62-135.69	0.024-0.040	135.95-136.00	0.033-0.046	135.90-135.93	0.023-0.036
16	139.62-139.69	0.031-0.037	139.95-140.01	0.001-0.041	139.92-139.94	0.023-0.038
17	143.71-143.78	0.018-0.043	144.08-144.14	0.010-0.044	144.04-144.08	0.029-0.040
18	147.81-147.88	0.028-0.039	148.19-148.24	0.020-0.042	148.14-148.18	0.033-0.039
19	151.93-152.01	0.024-0.038	152.31–152.37	0.018-0.048	152.27-152.30	0.031-0.042
D19S43	3					
6	172.80-172.89	0.025-0.039	173.29-173.39	0.013-0.043	173.28-173.33	0.028-0.034
7	176.77-176.85	0.030-0.038	177.25-177.35	0.003-0.047	177.25-177.30	0.035-0.042
8	180.72-180.82	0.028-0.042	181.25-181.35	0.004-0.047	181.25-181.30	0.023-0.040
9	184.79-184.92	0.019-0.043	185.37-185.47	0.013-0.045	185.37-185.44	0.031-0.040
10	188.80-188.92	0.028-0.043	189.42-189.51	0.021-0.042	189.40-189.47	0.035-0.042
11	192.70-192.81	0.023-0.047	193.33-193.42	0.010-0.053	193.31-193.39	0.028-0.045
12	196.67-196.79	0.033-0.044	197.32-197.44	0.007-0.045	197.32-197.40	0.029-0.046
12.2	198.81-198.94	0.025-0.046	199.44-199.56	0.001-0.046	199.46-199.53	0.038-0.048
13	200.64-200.75	0.035-0.043	201.28-201.40	0.005-0.043	201.28-201.36	0.028-0.048
13.2	202.61-202.71	0.030-0.045	203.26-203.36	0.010-0.045	203.26-203.33	0.027-0.043
14	204.53-204.64	0.033-0.045	205.19-205.29	0.016-0.052	205.19-205.26	0.037-0.040
14.2	206.51-206.62	0.023-0.047	207.17-207.27	0.019-0.052	207.16-207.23	0.029-0.043
15	208.44-208.54	0.031-0.048	209.11-209.21	0.008-0.050	209.10-209.18	0.028-0.046
15.2	210.41-210.50	0.029-0.050	211.08-211.19	0.009-0.050	211.09-211.16	0.020-0.044
16	212.36-212.45	0.027-0.051	213.03-213.13	0.003-0.048	213.03-213.11	0.037-0.039
16.2	214.34-214.43	0.032-0.047	215.03-215.14	0.031-0.048	215.03-215.11	0.032-0.045
17	216.42-216.52	0.033-0.048	217.12-217.25	0.023-0.046	217.15-217.23	0.029-0.044
17.2	218.34-218.44	0.027-0.049	219.05-219.19	0.003-0.062	219.07-219.15	0.035-0.048
18.2	222.38-222.48	0.031-0.044	223.12-223.25	0.017-0.054	223.14-223.22	0.036-0.055
19.2	226.31-226.42	0.036-0.057	227.06-227.22	0.015-0.061	227.11-227.19	0.034-0.052

	3130 <i>xl</i>		35	00	3500xL	
Allele	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
D22S10	145					
7	245.00-245.10	0.024-0.044	245.61-245.70	0.003-0.051	245.59-245.66	0.031-0.041
8	248.05-248.14	0.026-0.041	248.65-248.74	0.007-0.053	248.63-248.69	0.039-0.045
9	251.04-251.15	0.025-0.042	251.63-251.71	0.035-0.046	251.61-251.67	0.022-0.034
10	253.99-254.08	0.021-0.035	254.56-254.66	0.014-0.044	254.55-254.61	0.032-0.040
11	256.91-257.00	0.031-0.047	257.50-257.58	0.001-0.043	257.47-257.53	0.030-0.035
12	259.81-259.90	0.034-0.046	260.43-260.52	0.031-0.050	260.40-260.47	0.039-0.044
13	262.79-262.89	0.031-0.057	263.39-263.51	0.026-0.046	263.38-263.45	0.031-0.039
14	265.76-265.87	0.031-0.059	266.41-266.49	0.013-0.040	266.36-266.44	0.030-0.049
15	268.74-268.85	0.032-0.044	269.36-269.47	0.020-0.059	269.35-269.42	0.027-0.034
16	271.74-271.85	0.029-0.040	272.36-272.46	0.030-0.046	272.33-272.41	0.032-0.037
17	274.73-274.83	0.028-0.048	275.35-275.45	0.012-0.056	275.33-275.40	0.033-0.042
18	277.73-277.83	0.026-0.043	278.36-278.47	0.030-0.050	278.33-278.41	0.034-0.043
19	280.71-280.81	0.015-0.049	281.34-281.45	0.029-0.049	281.31-281.39	0.033-0.042
20	283.70-283.78	0.027-0.037	284.32-284.42	0.001-0.048	284.28-284.37	0.026-0.048
D10S12	248					
8	303.34-303.37	0.021-0.039	303.46-303.49	0.012-0.049	303.42-303.45	0.031-0.040
9	307.25-307.29	0.021-0.036	307.39-307.43	0.010-0.049	307.33-307.37	0.031-0.042
10	311.24-311.27	0.027-0.039	311.37-311.40	0.005-0.046	311.30-311.35	0.030-0.045
11	315.30-315.33	0.024-0.042	315.42-315.46	0.024-0.050	315.38-315.41	0.034-0.048
12	319.51-319.53	0.022-0.040	319.60-319.64	0.035-0.053	319.58-319.6	0.019-0.044
13	323.62-323.63	0.021-0.042	323.67-323.72	0.007-0.064	323.65-323.69	0.030-0.045
14	327.67-327.68	0.028-0.039	327.73-327.76	0.016-0.053	327.70-327.73	0.033-0.042
15	331.70-331.71	0.024-0.037	331.74-331.80	0.023-0.048	331.72-331.76	0.030-0.043
16	335.70-335.72	0.022-0.039	335.76-335.79	0.012-0.049	335.74-335.76	0.030-0.046
17	339.68-339.69	0.019-0.032	339.73-339.78	0.041-0.053	339.71-339.73	0.031-0.038
18	343.71-343.74	0.023-0.038	343.76-343.8	0.011-0.048	343.74-343.77	0.030-0.036
19	347.76-347.78	0.012-0.035	347.80-347.85	0.014-0.049	347.78-347.81	0.032-0.038

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 DNA Advisory Board (DAB) Standard 8.1.2.2).

Extra peaks: Stutter

Stutter definition

Stutter is a well-characterized PCR artifact that refers to the appearance of a minor peak one repeat unit smaller than the target STR allele product (minus stutter), or less frequently, one repeat larger (plus stutter) (Butler, 2005; Mulero *et al.*, 2006). Sequence analysis of stutter products at tetranucleotide STR loci has revealed that the minus stutter product is missing a single tetranucleotide core repeat unit relative to the main allele (Walsh *et al.*, 1996). Although 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.

Contact HID Support for more information on plus stutter.

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.

Stutter observations

Plus-stutter was regularly observed and was more significant in trinucleotide repeatcontaining loci (Figure 25 on page 83).

Peak heights were measured for amplified samples at the loci that are used in the kit. All data were generated on the 3500xL Genetic Analyzer. Some conclusions from these measurements and observations are:

- For each locus, the stutter percentage generally increases with allele length.
- Each allele within a locus displays a relatively consistent average stutter percentage.
- Peaks in the stutter position that are above the stutter filter percentage specified in the software are not filtered.
- The measurement of stutter percentage for allele peaks that are off-scale may be unusually high due to artificial truncation of the main allele peak.
- Stutter can be elevated when minus stutter and plus stutter overlap. This is typically observed when a given allele flanks another allele that is 2 repeat units away (as seen with the FGA locus in control 007 DNA).
- The magnitude and/or variability of stutter may increase with low DNA input amounts.

Figure 20 through Figure 24 show the stutter observed in the population study that are one repeat unit away from the alleles recorded. All data were generated on the 3500xL Genetic Analyzer.

The stutter filter settings that are derived from this data are listed in "Stutter percentage filter settings provided with GeneMapper" ID-X Software" on page 83.

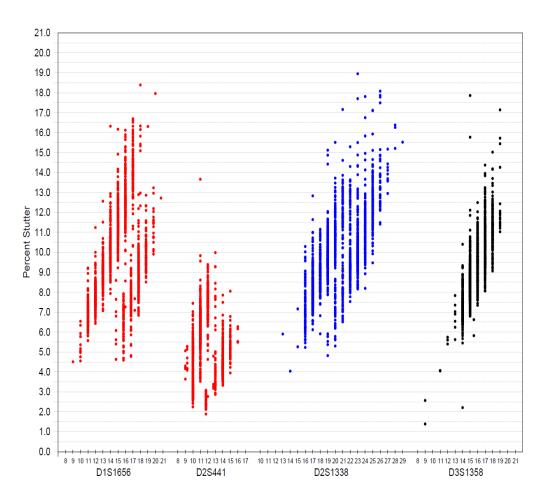


Figure 20 Stutter percentages for D1S1656, D2S441, D2S1338, and D3S1358 loci (Blue=FAM[™] dye, black=TED[™] dye, red=TAZ[™] dye)

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

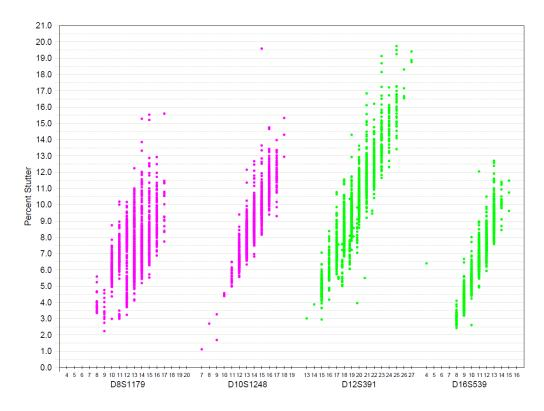


Figure 21 Stutter percentages for D8S1179, D10S1248, D12S391, and D16S539 (Green=VIC[™] dye and purple=SID[™] dye)

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

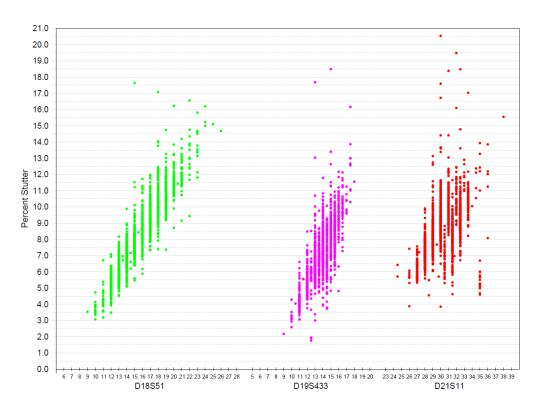


Figure 22 Stutter percentages for D18S51, D19S433, and D21S11 loci (Green=VIC[™] dye, red=TAZ[™] dye, purple=SID[™] dye)

 $\textbf{Note:} \ \ \text{The data in this figure are from the original formulation of the NGM Detect}^{\text{\tiny{M}}} \ kit.$

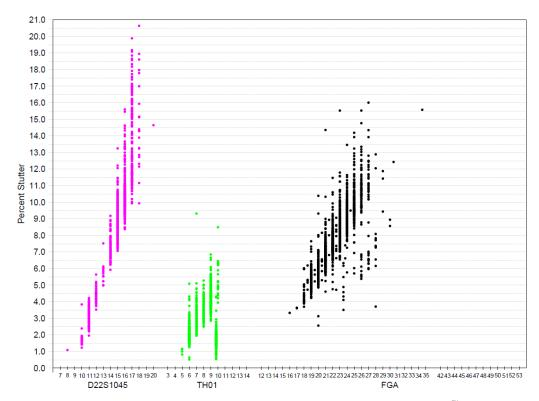


Figure 23 Stutter percentages for D22S1045, TH01, and FGA loci (Purple=SID[™] dye, Green=VIC[™] dye, black=TED[™] dye)

Note: Except for the TH01 marker data, the data in this figure are from the original formulation of the NGM $Detect^{\mathsf{TM}}$ kit.

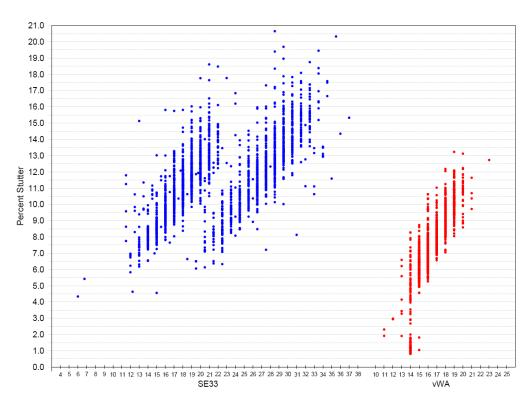


Figure 24 Stutter percentages for SE33 and vWA loci (Blue=FAM[™] dye, red=TAZ[™] dye)

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

Non-standard stutter peaks at the D22S1045 and SE33 loci

The D22S1045 locus in the NGM Detect[™] kit is a trinucleotide repeat locus, and shows an elevated level of plus stutter (Figure 25). Other loci, such as THO1 and D2S1338, may also exhibit relatively elevated plus stutter.

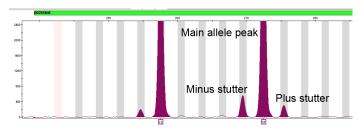


Figure 25 NGM Detect[™] kit electropherogram showing minus and plus stutter associated with the D22S1045 STR locus. Data produced on a 3500xL Genetic Analyzer.

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

STR loci such as D1S1656 and SE33 (Figure 26) contain more complex nucleotide sequences including regions of dinucleotide repeats which can yield additional stutter peaks. If these stutter peaks exceed the peak amplitude threshold (typically 175 RFU), they may be detected as additional alleles in the profile. The stutter file that is provided with the GeneMapper^{\mathbb{T}} *ID-X* Software for analysis of NGM Detect^{\mathbb{T}} kit data contains a minus 2–nt stutter filter for SE33 and D1S1656, as well as filters for commonly observed plus stutter, to prevent these peaks from being called in normal profiles.

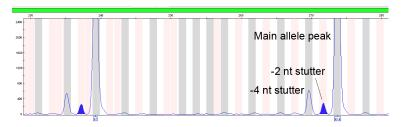


Figure 26 Example of a –2-nt reproducible stutter artifact at the SE33 locus. Data produced on a 3500xL Genetic Analyzer.

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

Stutter percentage filter settings provided with GeneMapper ${}^{\text{\tiny{M}}}$ /D-XSoftware

The settings in Table 5 and Table 6 were derived using the data that is shown in "Stutter observations" on page 77. The proportion of the stutter product relative to the main allele (stutter percent) is measured by dividing the height of the stutter peak by the height of the main allele peak.

Analysis showed that observed stutter data points were not normally distributed. As such, at each locus a best-fit, non-parametric statistical model was applied to the data and a threshold filter level that emulated historical stutter filter levels (approximately 99.7%) was derived.

IMPORTANT! The values that are shown in the table are the values that were determined during developmental validation studies using specific data sets. To determine the appropriate values to use for your applications, always perform internal validation studies.

Table 5 Minus stutter percentage filter settings provided with the GeneMapper[™] /D-X Software

Note: The data in this table are from the original validation study of the NGM Detect $^{\mathsf{TM}}$ kit. Based on the validation study with the updated kit formulation, no changes to the stutter filter values were required.

Locus ^[1]	% Stutter
D2S1338	17.61
SE33	18.63
SE33 (-2 nt)	6.41
D16S539	11.97
D18S51	16.15
TH01	7.66
D12S391	18.99
D3S1358	15.42
FGA	14.70
vWA	12.51
D21S11	16.88
D1S1656	16.68
D1S1656 (-2 nt)	3.90
D2S441	9.55
D8S1179	13.80
D19S433	13.57
D22S1045	19.02
D10S1248	14.32

^[1] These percentages are used as stutter filters in NGM_Detect_Stutter.txt

Table 6 Plus stutter percentage filter settings provided with the GeneMapper [™] *ID-X* Software

Note: The data in this table are from the original validation study of the NGM Detect $^{\text{TM}}$ kit. Based on the validation study with the updated kit formulation, no changes to the stutter filter values were required.

Locus ^[1]	% Stutter
D2S1338	11.55
SE33	9.80
D16S539	4.68
D18S51	8.19
TH01	6.14
D12S391	7.76
D3S1358	4.36
FGA	9.16
vWA	7.59
D21S11	9.14
D1S1656	5.47
D2S441	4.92
D8S1179	5.39
D19S433	6.91
D22S1045	8.22
D10S1248	2.60

^[1] These percentages are used as stutter filters in NGM_Detect_Stutter.txt

Extra peaks: Addition of 3' A nucleotide

3' A nucleotide addition definition

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 nontemplate 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.

3' A observations

The efficiency of +A addition is related to the particular sequence of the DNA at the 3′ end of the PCR product.

The NGM Detect[™] kit includes two main design features that promote maximum +A addition:

- The primer sequences have been optimized to encourage +A addition.
- The PCR chemistry allows complete +A addition with a short final incubation at 60°C for 5 minutes.

This final extension step gives the DNA polymerase additional time to complete +A addition to all double-stranded PCR products. Figure 27 shows examples of incomplete and normal +A addition. Final extension incubation for longer than the recommended time can result in double +A addition, in which two nontemplate adenosine residues are added to the PCR product. Double +A addition can cause "shoulders" on the right side of main allele peaks.

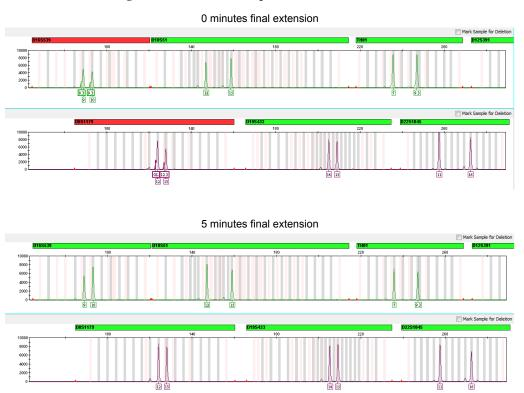


Figure 27 Omitting the final extension step results in shoulders on main allele peaks due to incomplete +A nucleotide addition. Examples shown are the smaller amplicons of VIC^{∞} and SID^{∞} dye channel data from a 3500xL Genetic Analyzer using the NGM Detect kit.

If the amount of input DNA is greater than recommended concentration, "shouldering" of allele peaks can be observed. Amplification of excess input DNA can also result in off-scale data and lowered IQCL peak heights. In this situation, the IQCL may also exhibit some shouldering due to incomplete +A nucleotide addition.

Extra peaks: Artifacts

Artifact definition

Artifacts and anomalies are seen in all molecular biological systems. Artifacts are typically reproducible. Anomalies are non-reproducible, intermittent occurrences that are not observed consistently in a system (for example, spikes and baseline noise).

Dye artifact observation

Due to improvements in PCR primer manufacturing processes, the incidence of artifacts has been greatly reduced in the NGM Detect[™] kit. Internal population studies show that kit electropherograms are free of reproducible dye artifacts in the kit read region of 64–458 nt. Two exceptions are as follows:

- A low level 113–117 nt dye artifact in the VIC[™] dye channel that has been detected below commonly used analytical thresholds.
- A low level ~66 nt dye artifact in the TED[™] dye channel. This artifact was approximately 40–80 RFU in our studies. The peak height observed may vary depending on the sensitivity of individual CE instruments.

Figure 28 shows the low baseline-level fluorescence that is observed in a typical negative control PCR. However, it is important to consider noise and other amplification-related artifacts when interpreting data.

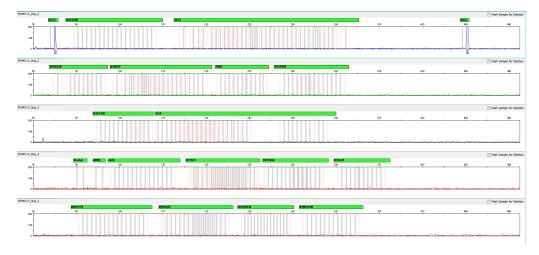


Figure 28 Examples of fluorescence background in data produced on a 3500xL Genetic Analyzer (Y-axis scale 0 to 200 RFU).

Some small PCR artifacts were occasionally observed with the human male 007 DNA provided as the control DNA in the NGM Detect $^{\text{\tiny M}}$ kit. The most prominent and consistent artifacts were:

- A 170-bp artifact in the SID[™] channel
- Two artifacts consistent with the size of a –2 and –3 repeat unit stutter for allele 14 of D2S441

In all cases, the peak heights of the artifacts were typically 1% or less than peak heights of the nearest true allele. If higher amounts of DNA (that is, 1 ng or more) are amplified, or if a lower PAT is used, then the peak heights could theoretically exceed the analytical threshold. The peak height of the artifacts appeared to be proportional to the amount of input 007 DNA in the PCR.

Characterization of loci

SWGDAM guideline 3.1

"The basic characteristics of a genetic marker should be determined and documented." (SWGDAM, December 2016)

Loci in this kit

This section describes basic characteristics of the 16 autosomal STR loci, Y indel locus, and sex-determining marker (Amelogenin), that are amplified with the NGM Detect $^{\text{\tiny M}}$ kit. Most of these loci have been extensively characterized by other laboratories.

Nature of polymorphisms

The primers for the Amelogenin locus flank a 6-nucleotide deletion in intron 1 of the X homolog. Amplification generates 108-nt and 114-nt products from the X and Y chromosomes, respectively. The primers for the Y indel flank a region in the q arm of the Y chromosome (Yq11.221). Depending on the haplotype of the sample, the amplification generates either a 91-nt or a 96-nt product. (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.) Except for D22S1045, a trinucleotide STR, the remaining loci are tetranucleotide short tandem repeat (STR) loci. The length differences among alleles of a particular locus are caused by differences in the number of repeat units.

We have sequenced all the alleles in the NGM Detect[™] kit Allelic Ladder, including microvariants. 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 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 Detect[™] 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; and Lareu, *et al.*, 1996).

Genetic linkage

Two sets of STR loci in the NGM Detect[™] 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; 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 1,034 individuals of three ethnic groups (350 African American, 349 Caucasian, and 335 Hispanic). 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 7 for results.

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

An independent analysis of data from the same collection of population samples (Budowle, *et al.*, 2010) also concluded that the 15 STR loci that are shared between the NGM[™] and NGM SElect[™] kits were independent at the population level (note that the SE33 locus was not part of this analysis). Therefore, to calculate the rarity of a profile for comparison to single-source and mixture samples, the frequencies of all loci including vWA and D12S391 could be multiplied. However, the analysis of the CEPH pedigree families demonstrated a degree of linkage between vWA and D12S391 that does not support the assumption of independence for kinship analysis.

Locus	Chromosome map position ^[1]	Chromosome Nuclear Coordinates ^[1] (million bp)	African- American (n = 350)	Caucasian (n = 350)	Hispanic (n = 293)
vWA	12p13.31	5.9	0.07	0.20	0.07
D12S391	12p13.2	12.2	0.86	0.29	0.27
D2S441	2p14	68	0.11	0.22	0.10
D2S1338	2q35	218	0.11 0.32	0.32	0.19

 Table 7
 GenePop software LD Result (p-value for pairwise analysis of loci)

Species specificity

SWGDAM Guideline 3.2

"The ability to detect genetic information from non-targeted species (e.g., detection of microbial DNA in a human assay) should be determined. The detection of genetic information from non-targeted species does not necessarily invalidate the use of the assay, but may help define the limits of the assay." (SWGDAM, December 2016)

Nonhuman studies

The NGM Detect $^{\text{m}}$ kit provides the required specificity for detecting human alleles. Species specificity testing was performed to ensure that there is no cross-reactivity with nonhuman DNA that may be present in forensic casework samples.

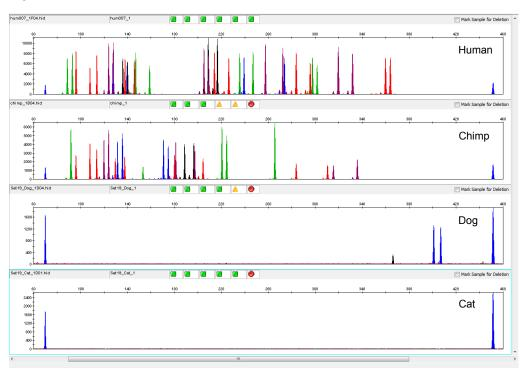
The following species were tested (in the specified amounts) using standard PCR and capillary electrophoresis conditions for the NGM DetectTM kit kit:

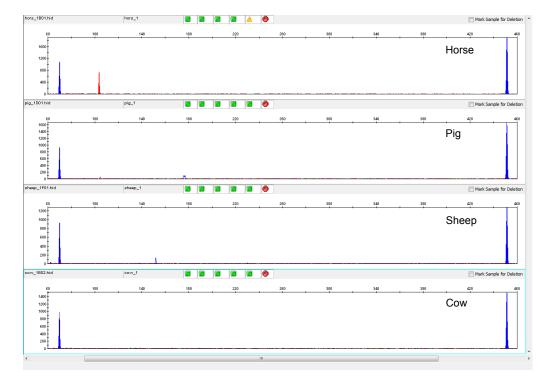
- Primates: gorilla, chimpanzee, orangutan, and cynomolgous (macaque) (0.5 ng each)
- Non-primates: mouse, dog, sheep, pig, rabbit, cat, horse, hamster, rat, chicken, and cow (5.0 ng each)
- Microorganisms: Candida albicans, Staphylococcus aureus, Escherichia coli, Neisseria gonorrhoeae, Bacillus subtilis, and Lactobacillus rhamnosus (pooled genomic DNAs, with approximately 50,000 copies of DNA from each species, per reaction)

^[1] 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.

Results were evaluated for the presence of any amplified peaks that would indicate cross reactivity of the NGM $Detect^{TM}$ kit with any of these non-human species.

Figure 29 shows select electropherogram results from the species specificity tests. The chimpanzee, gorilla (data not shown), orangutan (data not shown), and macaque (data not shown) DNA samples produced partial profiles in the 70–400 nucleotide region.





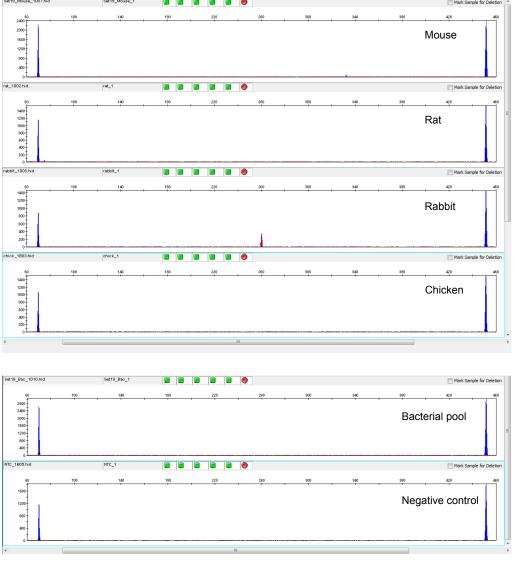


Figure 29 Representative electropherograms for some species tested in a species specificity study. Data produced on a 3500xL Genetic Analyzer.

Table 8 shows the most significant cross-reactive peaks that were observed among non-human, non-primate, genomic DNAs (that is, peaks over a 175 RFU Peak Amplitude Threshold on the 3500xL Genetic Analyzer). Peaks were observed for dog, horse, and rabbit. Most peaks did not fall into human STR locus bins or marker ranges, and would therefore not be confused with human STR alleles. This data shows that the likelihood of obtaining an allelic profile consistent with that from a human sample, from non-primates or microorganisms, is low.

Table 8 Observed cross-reactive peaks for non-human, non-primate animals.

Species	Dye channel	Size	Peak height
	D	401 bp	1,294 RFU
Dog	В	407 bp	1,046 RFU
	Υ	366 bp	289 RFU
Horse	R	104 bp	915 RFU
Rabbit	Р	260 bp	389 RFU

Sensitivity

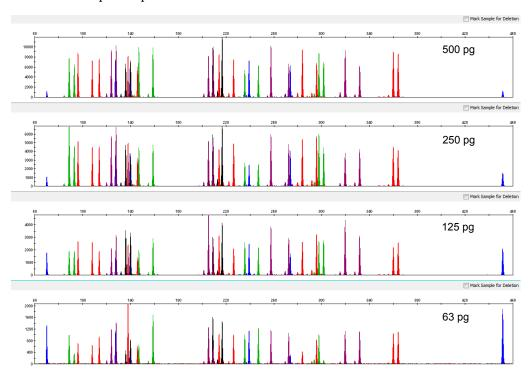
SWGDAM guideline 3.3

"The ability to obtain reliable results from a range of DNA quantities, to include the upper and lower limits of the assay, should be evaluated." (SWGDAM, December 2016)

Sensitivity observation

The recommended amount of input DNA for the NGM Detect^T kit is 0.5 ng for 30 cycles of amplification based on real-time PCR quantification, such as with the Quantifiler^T Trio DNA Quantification Kit or the Quantifiler HP DNA Quantification Kit. To determine the optimum input DNA, perform studies according to the quantification kit that you use.

If the sample contains degraded or inhibited DNA, amplification of a higher amount of DNA may be beneficial. In Figure 30, DNA Control 007 was serially diluted from 0.5–0.008 ng. Full profiles (35 PCR products) were consistently obtained at 0.125 ng, but occasional partial profiles resulted at lower concentrations.



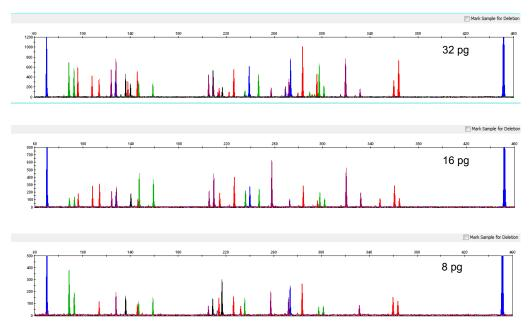


Figure 30 Electropherograms for amplifications using 500, 250, 125, 63, 32, 16, and 8 pg of DNA Control 007. Electrophoresis was performed on a 3500xL Genetic Analyzer. Note that as the DNA input is serially diluted by 2-fold, the Y-axis scale is also adjusted, as needed, to accommodate lower peak heights.

Stability

SWGDAM guideline 3.4

"The ability to obtain results from DNA recovered from biological samples deposited on various substrates and subjected to various environmental and chemical insults should be evaluated. 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 to determine the effects of such factors." (SWGDAM, December 2016)

Degraded DNA

As the average size of degraded DNA approaches the size of the target sequence, the amount of PCR product that is generated is reduced, because of the reduced number of intact templates in the size range required for amplification. Degraded DNA was prepared to examine the potential for differential amplification of loci. Approximately 10 μ g of high molecular weight DNA was sonicated and subjected to digestion with 0.25 U/ μ L of DNase I for various lengths of time (from ~5–10 minutes) to produce a range (low, medium, and high) of degradation levels (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 0.5 ng of degraded DNA using the NGM Detect 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.

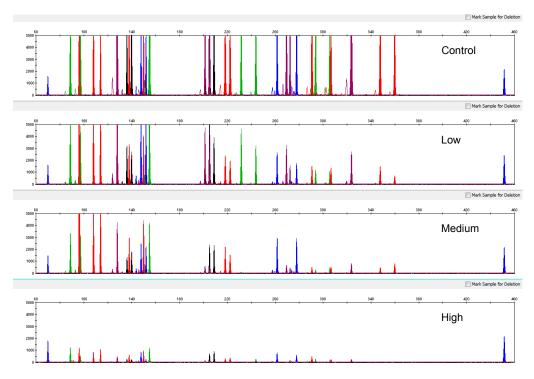


Figure 31 Amplification of a single donor DNA sample sonicated and incubated with $0.25 \text{ U/}\mu\text{L}$ of DNase I for various lengths of time (from ~5–10 minutes).

Effect of inhibitors: hematin

Heme compounds have been identified as PCR inhibitors in DNA samples that are extracted from bloodstains (DeFranchis *et al.*, 1988; Alkane *et al.*, 1994). It is believed that the inhibitor is co-extracted and co-purified with the DNA and then interferes with PCR by inhibiting polymerase activity. To examine the effects of hematin on the performance of the NGM DetectTM kit, 0.5 ng of DNA Control 007 was amplified in the presence of increasing concentrations of hematin for 30 cycles of amplification (Figure 32). The final concentrations of hematin that is used in the reaction were 0, 450, 575, and 700 μ M (Table 9).

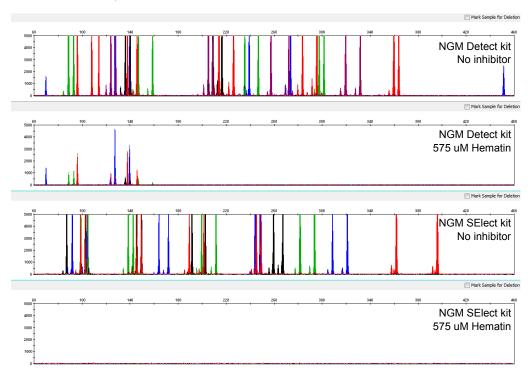


Figure 32 Electropherograms for the NGM Detect[™] kit and the NGM SElect[™] kit show the improved performance of the NGM Detect[™] kit in the presence of hematin compared with previous $AmpF\ell STR^{™}$ kits. The same set of inhibited samples was analyzed with the NGM Detect[™] kit and the NGM SElect[™] kit for 30 cycles of amplification.

Table 9 NGM Detect[™] kit performance in simulated hematin inhibition (N = 5).

Hematin concentration	Average number of alleles detected ^[1]	Minimum, maximum
0 μΜ	35	35, 35
450 μM	35	35, 35
575 μΜ	12.4	12, 13
700 μM	0	0, 0

^[1] Only those peaks ≥175 RFU were counted. A complete profile with DNA Control 007 yields 35 peaks using the NGM Detect[™] kit.

Effect of inhibitors: humic acid

Traces of humic acid may inhibit the PCR amplification of DNA evidence that is collected from soil. Amplification of 0.5 ng of DNA Control 007 in the presence of increasing amounts of humic acid was performed using the NGM Detect kit for 30 cycles of amplification (Figure 33). The final concentrations of humic acid tested in the reactions were 0, 140, 250 and 400 ng/ μ L (Table 10).

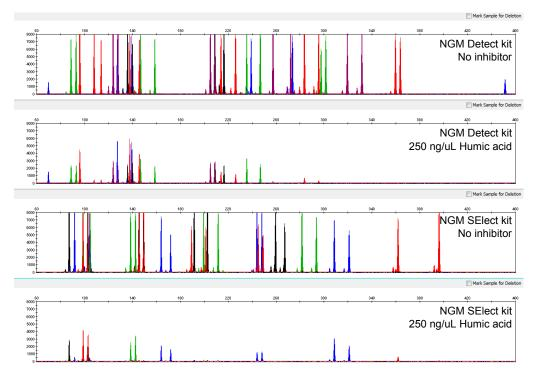


Figure 33 Electropherograms for the NGM Detect[™] kit and the NGM SElect[™] kit show the improved performance of the NGM Detect[™] kit in the presence of humic acid compared with previous $AmpF\ell STR^{™}$ kits. The same set of inhibited samples was analyzed with the NGM Detect[™] kit and the NGM SElect[™] kit for 30 cycles of amplification.

Table 10 NGM Detect[™] kit performance in simulated humic inhibition (N = 5).

Humic acid concentration	Number of alleles detected ^[1]	Minimum, maximum
0 ng/μL	35	35, 35
140 ng/μL	35	35, 35
250 ng/μL	26	26, 26
400 ng/μL	1	1, 1

^[1] Only those peaks ≥175 RFU were counted. A complete profile with DNA Control 007 yields 35 peaks using the NGM Detect[™] kit.

Mixture studies

SWGDAM guideline 3.8

"The ability to obtain reliable results from mixed source samples should be determined." (SWGDAM, December 2016)

Mixture study overview

Evidence samples may contain DNA from more than one individual. The possibility of multiple contributors should be considered when interpreting the results. Perform studies to determine a minimum peak height threshold to avoid typing when stochastic effects are likely to interfere with accurate interpretation of mixtures.

Evidence samples that contain body fluids and/or tissues originating from more than one individual are an important category of forensic casework.

It is essential to ensure that the DNA typing system is able to detect DNA mixtures. Typically, mixed samples can be distinguished from single-source samples by:

- The presence of more than two alleles at one or more loci
- 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.

If an unusually low peak height ratio is observed for one locus, and there are no other indications that the sample is a mixture, reamplify and reanalyze the sample 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

Mixture study observation

Median, minimum, and maximum peak height ratios observed for alleles in the NGM DetectTM kit loci in unmixed human population database samples (N=60) are shown in Figure 34. The donor samples that were used are listed in "Population samples used in these studies" on page 102.

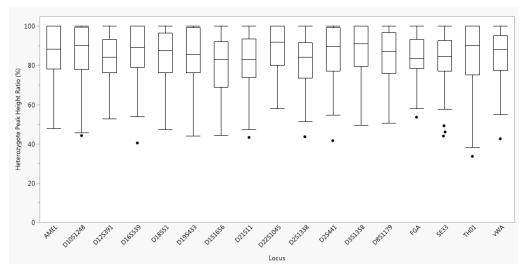


Figure 34 Heterozygote ratios for 0.5 ng of input DNA. The distribution of intra-locus peak height ratio is expressed as a percentage, by locus (low peak height allele divided by high peak height allele). Boxes show the middle 50% or interquartile range (IQR). Box halves below and above median show the second and third quartile, respectively. "Whiskers" indicate 1.5 IQR from the upper and lower margins of the IQR. Black dots are outlier data points more than 1.5 IQR from the median (N=60).

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 components at a single locus.

The ability to obtain and compare quantitative values for the different allele peak heights on Applied Biosystems^m 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 samples.

Note: Peak detection is a complex process that involves the STR chemistry, capillary electrophoresis conditions, and the data analysis software. Contact HID Support for a Technical Note with additional information on detecting peaks in electropherograms.

Limit of detection of the minor component

Mixtures of two DNA samples were examined at various ratios (0:1, 1:1, 1:3, 1:7, 1:15, 1:30, 30:1, 15:1, 7:1, 3:1 and 1:0 M:F ratios). The total amount of genomic input DNA mixed at each ratio was 0.5 ng. The samples were amplified in a GeneAmp^{$^{\text{TM}}$} PCR System 9700, then electrophoresed and detected using a 3500xL Genetic Analyzer.

The results of the mixed DNA samples are shown in Figure 35. The male 007 and female 9947a 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 7:1 (0.438 ng of 007 and 0.063 ng of 9947a) and 1:7 (0.063 ng of 007 and 0.438 ng of 9947a, with 2/3 replicates giving full profiles of the 007 minor contributor). 15:1, 30:1, 1:30, and 1:15 ratios resulted in partial profiles for the minor component. The genotypes of these samples are shown in Table 11.

Table 12 shows mixture sample compositions and STR allele counts from NGM Detect[™] assays.

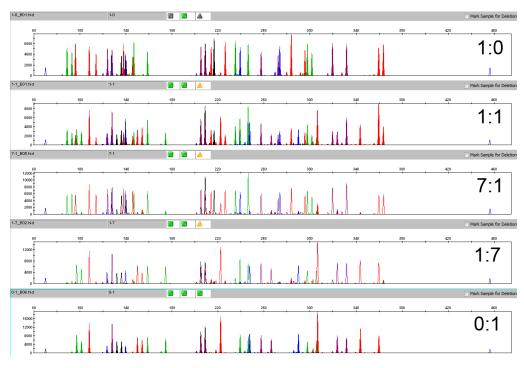


Figure 35 Amplification of DNA mixtures at various ratios. Panels show electropherograms for 1:0 (male 007 DNA only), 1:1 mixture, 7:1 mixture, 1:7 mixture, and 0:1 (female 9947a DNA only).

Note: The data in this figure are from the original formulation of the NGM Detect[™] kit.

Table 11 Genotypes of mixed DNA samples

Note: The data in this table are from the original validation study of the NGM Detect [™] kit.

Locus	Male 007 alleles		Female 99	47a alleles
IQCS	_	_	_	_
D2S1338	20	23	19	23
SE33	17	25.2	19	29.2
IQCL	_	_		_
D16S539	9	10	11	12
D18S51	12	15	15	19
TH01	7	9.3	8	9.3
D12S391	18	19	18	20
D3S1358	15	16	14	15
FGA	24	26	23	24
Y indel	2	_	_	_
Amelogenin	Х	Υ	X	
vWA	14	16	17	18
D21S11	28	31	30	_
D1S1656	13	16	18.3	_
D2S441	14	15	10	14
D8S1179	12	13	13	_
D19S433	14	15	14	15
D22S1045	11	16	11	14
D10S1248	12	15	13	15

Table 12Mixture sample compositions and STR allele counts from NGM Detect $^{™}$ assays

Note: The data in this table are from the original validation study of the NGM Detect $^{\text{TM}}$ kit.

Ratio (M:F)		ng per reaction	Number of unique alleles in non-stutter positions per contributor ^[1]		
	Male 007	Female 9947a	Total DNA	Male 007	Female 9947a
1:0	0.500	0.00	0.5	16.00	0.00
1:1	0.250	0.250	0.5	16.00	13.00
1:3	0.125	0.375	0.5	16.00	13.00
1:7	0.063	0.438	0.5	15.67	13.00

Ratio (M:F)		ng per reaction	Number of unique alle n non-stutter positions contributor ^[1]		
	Male 007	Female 9947a	Total DNA	Male 007	Female 9947a
1:15	0.031	0.469	0.5	12.33	13.00
1:30	0.016	0.484	0.5	8.00	13.00
30:1	0.484	0.016	0.5	16.00	8.67
15:1	0.469	0.031	0.5	16.00	11.67
7:1	0.438	0.063	0.5	16.00	13.00
3:1	0.375	0.125	0.5	16.00	13.00
0:1	0.000	0.500	0.5	0.00	13.00

^[1] Mixture sample compositions and STR allele counts from NGM Detect[™] assays. Average number of unique, distinct alleles (that is, alleles neither present in the other contributor genotype nor located in a stutter position) obtained in 3 replicate reactions per mixture sample. Full profiles of unique alleles in 007 and 9947a DNAs were 16 and 13, respectively.

Population data

SWGDAM guideline 3.7

"The distribution of genetic markers in populations should be determined in relevant population groups." (SWGDAM, December 2016)

Population data overview

To interpret the significance of a match between genetically typed samples, you must know the population distribution of alleles at each locus in question. If the genotype of the relevant evidence sample is:

- Different from the genotype of the reference sample for a suspect, then the suspect is *excluded* as the donor of the biological evidence that was tested. An exclusion is independent of the frequency of the two genotypes in the population.
- The same as the genotype of the reference sample for a suspect, then the suspect is *included* as a possible source of the evidence sample.

The probability that another, unrelated individual would also match the evidence sample is estimated by the frequency of that genotype in the relevant populations.

Loci in the kit

The NGM DetectTM kit contains loci for which extensive population data are available. For additional information on the loci shared between many of the AmpF ℓ STRTM kits, see the population data and additional studies section of the $AmpF\ell STR^{TM} NGM$ $SElect^{TM} PCR$ Amplification Kit Kit User Guide (Pub. No. 4458841) and the $AmpF\ell STR^{TM}$ $Identifiler^{TM} Plus$ PCR $STR^{TM} Identifiler^{TM} Plus$ $STR^{TM} Identifiler^{TM} Plus$ S

Population samples used in these studies

Note: The data in this section are from the original validation study of the NGM Detect^{M} kit.

The NGM Detect[™] kit has high genotypic concordance (>99.5%) to the NGM SElect[™] kit and by extension to the GlobalFiler[™] kit. (All STR markers in the NGM SElect[™] kit are included in the GlobalFiler[™] kit with both the GlobalFiler[™] and NGM SElect[™] kits sharing the same primers for common loci.) Therefore, the existing population data from the GlobalFiler[™] kit was used, as applicable, to generate the population data that is provided in this section. Whole blood samples, provided by the Interstate Blood Bank (Memphis, Tennessee) and Boca Biolistics (Coconut Creek, Florida), were collected in the United States (with no geographical preference) from randomly selected individuals of known ethnicities. Ethnicities of sample donors were:

- African-American—330 samples
- Asian—153 samples
- Caucasian—343 samples
- Hispanic—368 samples

DNA was extracted using a 6100 Nucleic Acid Prep Station.

In addition to the alleles that we 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).

Concordance studies

In the validation study of the original NGM Detect[™] kit formulation, a concordance population study with 1,165 in-house samples identified two samples that were discordant because of allele dropout or severe allele imbalance in the NGM Detect assay. The samples were characterized and degenerate oligos were added to the SE33 (2 oligos) and D8S1179 (1 oligo) markers in the final stages of development. The resulting concordance between the NGM SElect kit and the NGM Detect kit was greater than 99.5%.

In the validation study of the updated NGM Detect[™] kit formulation, we used the new primer set with 1,092 individuals from the same donor population. We observed 100% concordance between the new primer set and the original primer set at the TH01 locus.

Probability of Identity definition

The P_I value is the probability that two individuals selected at random will have an identical genotype (Sensabaugh, 1982).

Probability of identity

Table 13 shows the Autosomal STR allele frequencies at NGM Detect $^{\text{\tiny TM}}$ kit loci by population group.

Table 14 shows the Y-specific allele frequencies by population group for NGM Detect^T kit Y indel locus. The Y-specific allele frequencies were not included in the probability of identity calculation.

Table 15 shows the Probability of identity (P_I) values of the NGM DetectTM kit loci individually and combined.

Table 13 Autosomal allele frequencies by population group for NGM Detect[™] kit STR loci. (*=Alleles not detected or not detected in significant quantities)

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)		
D2S1338						
10	*	*	*	*		
11	*	*	*	*		
12	*	*	*	*		
13	0.15*	*	0.15*	*		
14	*	*	0.15*	*		
15	0.30*	*	0.15*	*		
16	5.3	1.63	4.08	3.8		
17	10	14.05	18.37	17.8		
18	4.85	13.07	8.31	6.52		
19	16.21	16.67	14.14	17.53		
20	10.45	8.82	15.74	13.86		
21	11.97	2.94	2.92	3.67		
22	12.42	5.88	1.75	6.52		
23	9.24	18.3	10.06	14.27		
24	8.79	11.11	10.2	8.83		
25	6.97	5.88	12.1	5.43		
26	2.58	*	1.6	1.49		
27	0.76	0.33*	0.29*	0.14*		
28	*	0.98*	*	0.14*		
29	*	0.33*	*	*		

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
SE33				
3.2	*	*	*	*
4.2	*	*	*	*
5	*	*	*	*
6	*	*	*	*
6.3	*	*	0.15*	*
7	*	*	*	*
8	*	*	*	*
8.2	*	*	*	*
9	*	*	*	*
9.2	*	*	*	*
10	*	*	*	*
10.2	*	*	*	*
11	*	*	*	*
11.2	0.76	*	*	0.14*
12	0.15*	0.33*	0.44*	0.14*
12.1	*	*	0.15*	*
12.2	0.30*	*	0.15*	0.14*
13	1.36	*	0.87	1.22
13.2	0.45*	*	*	0.14*
14	3.33	*	3.64	1.77
14.2	0.15*	*	*	0.82
14.3	*	*	0.15*	*
15	4.24	1.31*	3.64	4.89
15.2	0.15*	*	0.15*	0.14*
15.3	*	*	*	*
16	6.97	3.59	5.39	5.57
16.2	0.30*	*	*	0.27*
16.3	*	*	*	0.14*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
17	7.73	5.23	6.56	8.7
17.2	0.15*	*	*	*
17.3	*	*	0.15*	*
18	10.76	4.9	7.87	10.05
18.2	0.15*	*	*	0.27*
19	15	9.48	8.31	8.15
19.2	0.30*	*	0.29*	*
19.3	*	*	0.29*	*
20	9.55	6.86	5.25	4.48
20.2	0.91	0.33*	0.87	0.82
21	5.76	6.21	2.04	3.12
21.1	*	*	*	*
21.2	0.91	1.63	1.17	1.09
22	1.97	2.61	0.58*	1.09
22.2	1.36	2.29	3.35	2.31
23	0.30*	*	*	*
23.2	0.61*	2.61	2.62	2.85
23.3	*	*	*	0.14*
24	0.30*	0.33*	0.15*	0.14*
24.2	1.67	6.54	4.52	2.31
25	*	*	*	*
25.2	2.42	7.19	3.79	3.12
26	0.15*	*	*	0.27*
26.2	5.61	7.52	4.52	6.39
27	*	*	*	*
27.2	5.91	3.59	6.85	7.07
28	*	*	*	*
28.2	3.94	7.84	7.73	6.25
29	*	*	*	*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
29.2	2.58	8.5	7.87	5.84
30	*	*	*	*
30.2	1.21	7.52	4.66	3.8
31	*	*	*	0.14*
31.2	1.06	1.63	2.77	2.31
32	*	*	0.44*	*
32.2	0.76	1.31*	1.6	2.04
33	*	*	0.29*	0.41*
33.2	0.45*	0.33*	0.15*	0.54*
34	*	*	0.29*	0.41*
34.2	0.15*	*	*	0.27*
35	*	*	0.15*	*
35.2	*	0.33*	*	*
36	*	*	0.15*	*
36.2	*	*	*	*
37	*	*	*	0.14*
37.2	*	*	*	*
38	*	*	*	*
38.2	*	*	*	*
39	*	*	*	*
39.2	*	*	*	*
40	*	*	*	*
41	*	*	*	*
42	*	*	*	*
43	*	*	*	*
D16S539		'	· 	
4	*	*	*	*
5	*	*	*	*
6	*	*	*	0.14*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
7	*	*	*	*
8	3.33	*	1.46	2.04
8.3	*	*	*	*
9	21.67	31.05	12.68	10.19
9.3	*	*	*	*
10	11.52	14.05	4.08	15.76
11	30	20.59	32.22	31.79
11.3	*	*	*	*
12	19.09	21.57	30.9	24.18
12.1	*	*	*	*
12.2	*	*	*	*
13	13.03	11.44	16.76	14.4
13.3	*	*	*	*
14	1.36	1.31*	1.75	1.22
15	*	*	0.15*	0.27*
16	*	*	*	*
D18S51				
6	*	*	*	*
7	*	*	*	*
8				
9	*	*	*	0.14*
9.2	*	*	*	*
10	0.15*	*	1.17	0.68
10.2	0.15*	*	*	*
11	0.45*	1.31*	0.87	1.22
11.2	*	*	*	*
12	6.21	5.56	15.01	10.46
12.2	*	*	*	*
12.3	*	*	*	*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
13	3.94	17.32	11.95	11.41
13.1	*	*	*	*
13.2	0.30*	*	*	*
13.3	*	*	*	*
14	5.91	22.88	17.64	16.3
14.2	0.45*	*	*	0.14*
15	16.52	16.99	15.31	12.23
15.2	*	*	*	0.14*
15.3	*	*	*	*
16	18.18	12.42	11.95	12.91
16.1	*	*	*	*
16.2	*	*	*	*
17	16.36	6.54	10.79	17.39
17.2	*	*	*	*
17.3	*	*	*	*
18	14.09	4.9	8.31	7.74
18.1	*	*	*	*
18.2	*	*	*	*
19	9.7	5.23	4.08	3.53
19.2	*	*	*	*
20	4.7	1.96	1.31	1.9
20.2	0.15*	*	*	*
21	1.82	1.96	1.02	2.17
21.2	*	*	*	*
22	0.61*	0.98*	0.29*	0.68
22.2	*	*	*	*
23	0.30*	0.98*	0.29*	0.54*
23.2	*	*	*	*
24	*	0.65*	*	0.27*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
25	*	*	*	0.14*
26	*	0.33*	*	*
27	*	*	*	*
28	*	*	*	*
TH01				
3	*	*	*	*
4	*	*	*	*
5	0.45*	*	0.15*	*
5.3	*	*	*	*
6	15.45	13.07	21.72	27.17
6.3	*	*	*	*
7	37.42	26.14	17.64	32.74
7.3	*	*	*	*
8	20.61	3.59	11.37	8.7
8.3	*	*	*	*
9	16.06	51.63	17.06	12.77
9.3	8.33	4.25	31.2	17.12
10	1.52	1.31*	0.87	1.49
10.3	*	*	*	*
11	*	*	*	*
12	*	*	*	*
13	*	*	*	*
13.3	*	*	*	*
14.3	*	*	*	*
D12S391				
13	*	*	*	0.14*
14	*	*	*	0.14*
15	7.58	2.61	4.37	4.08
16	5.15	0.98*	3.35	5.03

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
17	16.52	8.17	10.35	7.34
17.1	0.45*	*	*	0.27*
17.3	0.61*	*	1.9	1.22
18	24.55	28.43	16.18	19.7
18.1	*	*	*	*
18.3	1.21	*	2.19	2.17
19	13.94	23.86	12.54	18.75
19.1	0.61*	*	*	*
19.3	0.30*	*	0.58*	1.22
20	11.52	16.99	9.77	17.12
20.1	*	*	*	*
20.3	*	*	0.15*	*
21	7.27	10.78	13.56	8.7
21.3	0.15*	*	0.15*	*
22	5	3.59	10.79	6.79
23	3.64	3.27	8.16	3.67
24	0.61*	0.98*	3.64	1.9
24.3	*	*	*	*
25	0.61*	0.33*	1.9	1.36
26	*	*	0.29*	0.27*
27	*	*	0.15*	0.14*
D3S1358				
8	*	*	*	*
9	0.30*	*	*	0.14*
10	*	*	*	*
11	*	*	0.29*	*
12	0.15*	0.33*	*	0.14*
13	0.61*	*	0.15*	0.41*
14	9.09	2.61	15.16	9.1

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
15	28.18	49.02	27.26	34.65
15.2	0.30*	*	*	*
16	32.42	21.9	24.34	26.9
16.2	*	*	*	*
17	22.27	19.61	19.68	17.93
17.1	*	*	*	*
17.2	*	*	*	*
18	6.06	6.54	11.66	9.92
18.2	*	*	*	*
19	0.61*	*	1.46	0.82
20	*	*	*	*
20.1	*	*	*	*
21	*	*	*	*
FGA				
12.2	*	*	*	*
13	*	*	*	*
14	*	*	*	*
15	*	*	*	*
16	*	0.33*	0.15*	*
16.1	0.30*	*	*	*
16.2	*	*	*	*
17	*	0.33*	0.15*	*
17.2	*	*	*	*
18	0.91	3.27	1.02	0.68
18.2	0.61*	*	*	*
19	6.97	4.25	5.69	7.61
19.2	0.45*	*	*	*
19.3	*	*	*	*
20	6.82	3.92	14.87	8.7

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
20.1	*	*	*	*
20.2	0.30*	*	0.44*	0.27*
20.3	*	*	*	*
21	11.67	13.07	18.22	13.45
21.2	0.15*	0.33*	0.29*	*
21.3	*	*	*	*
22	17.27	14.38	19.24	14.4
22.1	*	*	*	*
22.2	0.15*	*	0.87	0.54*
22.3	*	*	*	*
23	17.27	27.12	14.87	12.91
23.1	*	*	*	*
23.2	*	0.65*	0.44*	0.41*
23.3	0.30*	*	*	*
24	18.94	18.3	14.43	15.62
24.1	*	*	*	*
24.2	*	0.33*	*	*
24.3	*	*	*	*
25	9.55	9.8	6.71	13.72
25.1	*	*	*	*
25.2	*	*	*	*
25.3	*	*	*	*
26	4.09	3.27	1.9	7.07
26.1	*	*	*	*
26.2	*	*	*	*
27	2.58	0.65*	0.58*	3.12
27.2	*	*	*	*
28	1.21	*	0.15*	0.95
28.2	*	*	*	*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
29	*	*	*	0.41*
29.2	*	*	*	*
30	0.15*	*	*	0.14*
30.2	0.15*	*	*	*
31	*	*	*	*
31.2	*	*	*	*
32	*	*	*	*
32.2	*	*	*	*
33.2	*	*	*	*
34.2	0.15*	*	*	*
41.2	*	*	*	*
42.2	*	*	*	*
43.2	*	*	*	*
44.2	*	*	*	*
45.2	*	*	*	*
46.2	*	*	*	*
47.2	*	*	*	*
48.2	*	*	*	*
49.2	*	*	*	*
50.2	*	*	*	*
51.2	*	*	*	*
52.2	*	*	*	*
vWA				
10	*	*	*	*
11	0.45*	*	*	0.14*
12	*	*	*	0.27*
13	0.91	*	0.15*	0.14*
14	7.27	23.53	8.75	6.52
15	20.91	1.63	12.24	9.78

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
15.2	*	*	*	*
16	27.58	15.36	22.3	30.57
17	19.85	29.74	27.41	27.17
18	13.79	19.61	17.78	18.07
18.2	*	*	*	*
19	6.52	9.15	10.06	6.39
20	1.97	0.98*	1.31	0.82
21	0.61*	*	*	*
22	*	*	*	*
23	0.15*	*	*	*
24	*	*	*	*
25	*	*	*	*
D21S11				
23	*	*	*	*
23.2	*	*	*	*
24	*	*	*	*
24.2	*	*	*	0.27*
24.3	*	*	*	*
25	*	*	*	*
25.2	*	*	*	*
25.3	*	*	*	*
26	0.30*	*	0.58*	0.41*
26.2	*	*	*	*
27	5.91	*	2.62	1.49
27.1	*	*	*	*
27.2	*	*	*	*
28	25.15	4.9	16.76	11.41
28.1	*	*	*	*
28.2	*	0.65*	*	0.14*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
28.3	*	*	*	*
29	15.61	26.8	23.76	21.06
29.1	*	*	*	*
29.2	*	*	0.15*	*
29.3	0.15*	*	0.15*	*
30	20.76	30.72	23.18	27.17
30.1	*	*	*	*
30.2	1.67	0.65*	2.77	1.77
30.3	*	*	*	*
31	8.79	9.48	6.85	5.16
31.1	*	*	*	*
31.2	4.55	3.92	8.89	11.14
31.3	*	0.33*	*	*
32	1.36	2.61	2.33	1.36
32.1	*	*	*	*
32.2	7.12	14.38	9.62	12.5
32.3	*	*	*	*
33	0.91	0.98*	*	0.14*
33.1	*	*	*	*
33.2	3.18	4.58	1.9	5.3
33.3	*	*	*	*
34	0.15*	*	*	*
34.1	*	*	*	*
34.2	*	*	0.44*	0.14*
35	3.64	*	*	0.27*
35.1	*	*	*	*
35.2	*	*	*	*
36	0.76	*	*	0.14*
36.1	*	*	*	*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
36.2	*	*	*	*
37	*	*	*	*
37.2	*	*	*	*
38	*	*	*	0.14*
38.2	*	*	*	*
39	*	*	*	*
D1S1656				
8	*	*	*	*
9	0.15*	*	*	0.14*
10	1.36	*	0.29*	0.41*
11	5.3	3.59	6.27	3.94
12	8.48	4.25	15.74	9.38
13	11.06	13.73	7	7.07
13.3	*	*	*	*
14	25	6.21	6.27	11.28
14.3	0.91	*	0.29*	0.27*
15	16.97	20.26	15.31	15.49
15.3	1.82	*	8.75	2.99
16	10	31.05	9.33	15.08
16.3	7.27	0.65*	4.96	5.16
17	2.73	14.05	4.96	6.79
17.1	*	*	0.29*	*
17.3	5.76	3.59	12.68	15.76
18	0.45*	0.33*	0.29*	0.82
18.3	1.82	1.63	5.98	4.48
19	0.15*	*	*	*
19.3	0.61*	0.33*	1.6	0.68
20	*	0.33*	*	*
20.3	0.15*	*	*	*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
21	*	*	*	*
21.3	*	*	*	*
D2S441				
7	*	*	*	*
8	0.15*	*	*	*
9	*	*	0.58*	0.14*
9.1	*	2.94	*	*
10	9.09	20.59	19.83	30.3
11	35.61	36.27	33.09	31.93
11.3	2.88	2.61	5.1	4.62
12	20.45	20.92	4.08	3.8
12.3	0.15*	*	0.29*	0.41*
13	3.48	6.21	3.35	1.9
13.3	*	*	*	*
14	26.21	9.8	28.86	23.1
14.3	*	*	*	*
15	1.97	0.65*	4.37	3.4
16	*	*	0.44*	0.41*
17	*	*	*	*
18	*	*	*	*
D8S1179				
4	*	*	*	*
5	*	*	*	*
6	*	*	*	*
7	*	*	*	*
8	0.30*	*	2.04	0.68
9	0.30*	*	1.31	0.27*
10	3.33	9.8	10.5	9.51
10.2	*	*	*	*

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
11	5.61	9.48	6.71	5.03
12	11.36	13.4	15.16	12.5
12.3	*	*	*	*
13	18.18	24.18	33.24	33.15
13.3	*	*	*	*
14	35.91	15.69	18.8	23.23
15	18.03	22.22	9.04	11.41
15.3	*	*	*	*
16	5.91	4.25	2.77	3.53
17	1.06	0.98*	0.44*	0.68
18	*	*	*	*
19	*	*	*	*
20	*	*	*	*
D19S433				
5	*	*	*	*
5.2	*	*	*	*
6	*	*	*	*
6.2	*	*	*	*
7	*	*	*	*
8	*	*	*	*
9	0.30*	0.33*	*	*
9.2	*	*	*	*
10	1.21	*	0.15*	0.41*
10.2	0.15*	*	*	*
11	9.85	*	*	1.63
11.2	0.30*	*	*	0.27*
12	10.45	4.58	7.29	8.42
12.1	*	*	0.15*	*
12.2	3.94	0.33*	0.15*	1.49

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
13	27.88	28.1	27.26	18.48
13.1	*	*	*	*
13.2	5.3	2.61	1.6	6.93
14	18.94	23.2	35.13	30.71
14.2	5.3	9.48	2.04	4.62
14.3	*	*	*	*
15	6.67	7.52	16.18	13.04
15.2	4.39	20.26	3.5	6.79
16	1.52	0.33*	5.69	4.08
16.2	3.18	2.61	0.29*	2.17
17	*	*	0.29*	0.54*
17.2	0.61*	0.65*	0.15*	0.41*
18	*	*	0.15*	*
18.2	*	*	*	*
19	*	*	*	*
19.2	*	*	*	*
20.2	*	*	*	*
D22S1045		1		
6	*	*	*	*
7	*	*	*	*
8	0.61*	*	*	*
9	*	*	*	*
10	4.09	*	0.44*	0.68
11	14.7	15.36	13.85	7.61
12	6.21	0.33*	0.58*	0.95
13	0.30*	0.33*	1.02	1.09
14	7.88	0.33*	3.35	2.04
15	23.33	33.66	36.3	43.48
16	20.3	23.86	36.3	34.65

Allele	African American (n=330)	Asian (n=153)	U.S. Caucasian (n=343)	U.S. Hispanic (n=368)
17	20.45	24.18	7.58	8.42
18	2.12	1.96	0.58*	0.95
19	*	*	*	*
20	*	*	*	0.14*
21	*	*	*	*
D10S1248				
7	0.15*	*	*	*
8	*	*	*	0.14*
9	0.15*	*	*	0.14*
10	*	0.33*	*	0.14*
11	3.64	*	0.58*	0.27*
12	14.09	10.78	3.5	4.48
13	22.88	36.93	29.45	25.95
14	27.88	22.55	29.74	36.14
15	18.48	22.55	19.39	22.69
16	10.15	5.23	13.41	7.74
17	2.27	1.63	3.64	2.31
18	0.30*	*	0.29*	*
19	*	*	*	*
20	*	*	*	*

Table 14 Y-specific frequencies by population group for the NGM Detect[™] kit Y indel locus. (*=Alleles not detected or not detected in significant quantities)

Allele	African American (n = 246)	Asian (n = 65)	U.S. Caucasian (n = 233)	U.S. Hispanic (n = 182)
Y indel				
1	1.22	67.69	*	0.55*
2	98.78	32.31	100	99.45

Table 15 Probability of identity (P_I) values for the NGM Detect[™] kit STR loci

Locus	African American (n = 330)	Asian (n = 153)	U.S. Caucasian (n = 343)	U.S. Hispanic (n = 368)
D10S1248	0.0693	0.1045	0.0943	0.1131
D12S391	0.0377	0.0664	0.0231	0.0318
D16S539	0.0727	0.0915	0.1043	0.0809
D18S51	0.0322	0.0402	0.0311	0.0281
D19S433	0.0388	0.0663	0.0862	0.0484
D1S1656	0.034	0.0564	0.0223	0.0247
D21S11	0.0453	0.0671	0.052	0.0487
D22S1045	0.0559	0.1073	0.1309	0.1604
D2S1338	0.0225	0.0337	0.0316	0.0316
D2S441	0.103	0.0961	0.0976	0.1079
D3S1358	0.0984	0.1689	0.0749	0.0949
D8S1179	0.0762	0.0527	0.0631	0.0661
FGA	0.0322	0.0555	0.0384	0.0282
SE33	0.0118	0.0139	0.0085	0.0081
TH01	0.0949	0.175	0.0801	0.0902
vWA	0.0622	0.084	0.065	0.0926
Combined	7.96 × 10 ⁻²²	2.25 × 10 ⁻¹⁹	2.29 × 10 ⁻²¹	3.12 × 10 ⁻²¹

Probability of paternity exclusion observation

Allele frequencies, observed heterozygosity (Ho), expected heterozygosity (He), Match Probability (MP), and p-value of each locus was calculated using a software program developed by Ge (Li *et al.*, 2013) and shown in the following table.

Departures from Hardy-Weinberg Equilibrium (HWE) expectations of each locus were derived using Arlequin (Excoffier *et al.*, 2010). After Bonferroni correction (Weir, 1990), 024), no departures from HWE were checked at any locus.

(p-value = 0.05/21 = 0.0024), no departures from HWE were observed at any locus.

Table 16 Allele frequencies, observed heterozygosity (Ho), expected heterozygosity (He), Match probability (MP), and p-value of STR loci

		African A	American	ı		As	ian			U.S. Ca	ucasian			U.S. H	spanic	
Marker	Но	He	MP	p- value	Но	He	MP	p- value	Но	Не	MP	p- value	Но	He	MP	p- value
Y-indel	-	-	0.976	-	-	-	0.556	-	-	-	1	-	-	-	1	-
D3S1358	0.772	0.762	0.094	0.902	0.686	0.682	0.151	0.012	0.753	0.786	0.079	0.698	0.698	0.758	0.098	0.291
vWA	0.752	0.798	0.068	0.116	0.829	0.784	0.08	0.309	0.841	0.807	0.064	0.621	0.842	0.782	0.08	0.463
D16S539	0.772	0.797	0.071	0.026	0.8	0.764	0.092	0.457	0.801	0.749	0.103	0.047	0.77	0.774	0.084	0.835
D8S1179	0.782	0.793	0.068	0.662	0.814	0.828	0.052	0.795	0.793	0.797	0.065	0.122	0.755	0.802	0.064	0.518
D21S11	0.861	0.849	0.039	0.553	0.8	0.791	0.069	0.667	0.873	0.837	0.046	0.385	0.827	0.839	0.044	0.315
D18S51	0.931	0.868	0.031	0.324	0.829	0.853	0.038	0.572	0.873	0.872	0.03	0.962	0.849	0.87	0.031	0.945
D2S441	0.772	0.756	0.099	0.421	0.714	0.746	0.101	0.238	0.757	0.766	0.09	0.077	0.791	0.763	0.094	0.611
D19S433	0.812	0.825	0.051	0.663	0.714	0.802	0.064	0.13	0.785	0.774	0.083	0.859	0.82	0.834	0.046	0.446
TH01	0.762	0.747	0.102	0.418	0.614	0.656	0.171	0.381	0.753	0.783	0.081	0.326	0.77	0.767	0.091	0.845
FGA	0.782	0.866	0.033	0.082	0.9	0.841	0.044	0.24	0.829	0.857	0.037	0.337	0.842	0.882	0.025	0.127
D22S1045	0.842	0.822	0.055	0.062	0.743	0.742	0.112	0.966	0.705	0.714	0.131	0.026	0.698	0.672	0.162	0.064
SE33	0.96	0.929	0.009	0.776	0.943	0.936	0.008	0.526	0.968	0.947	0.005	0.532	0.921	0.941	0.007	0.597
D10S1248	0.792	0.789	0.075	0.823	0.757	0.764	0.091	0.928	0.785	0.769	0.09	0.63	0.691	0.724	0.124	0.336
D1S1656	0.921	0.863	0.033	0.351	0.757	0.818	0.056	0.043	0.912	0.899	0.019	0.55	0.871	0.896	0.02	0.048
D12S391	0.861	0.864	0.032	0.19	0.771	0.808	0.063	0.65	0.904	0.896	0.02	0.45	0.842	0.874	0.028	0.071
D2S1338	0.911	0.894	0.02	0.763	0.871	0.872	0.03	0.356	0.88	0.878	0.027	0.23	0.906	0.877	0.027	0.929

The following table shows the Probability of paternity exclusion (PE) values of the NGM Detect $^{\text{\tiny TM}}$ kit STR loci individually and combined.

The PE value is the probability, averaged over all possible mother-child pairs, that a random alleged father will be excluded from paternity after DNA typing using the NGM Detect[™] kit STR loci (Chakraborty, Stivers, and Zhong, 1996).

Table 17 Probability of paternity exclusion values for the NGM Detect[™] kit STR loci

Locus	African American (n = 330)	Asian (n = 153)	Caucasian (n = 343)	Hispanic (n = 368)
D10S1248	0.6623	0.5353	0.5649	0.4644
D12S391	0.7401	0.631	0.8032	0.6588
D16S539	0.5548	0.6063	0.5915	0.5623
D18S51	0.7892	0.656	0.7557	0.7121
D19S433	0.6332	0.5238	0.5135	0.6431
D1S1656	0.7462	0.5703	0.8032	0.7338
D21S11	0.728	0.6063	0.7264	0.7013
D22S1045	0.7038	0.4795	0.4507	0.397
D2S1338	0.814	0.7463	0.7498	0.7392
D2S441	0.5228	0.5353	0.4986	0.5051
D3S1358	0.4918	0.3976	0.5338	0.4689
D8S1179	0.599	0.6063	0.6187	0.5381
FGA	0.728	0.8397	0.6632	0.7175
SE33	0.8639	0.88	0.9231	0.8781
TH01	0.5124	0.3424	0.5036	0.5381
vWA	0.6103	0.6186	0.6576	0.6276
PEi	7.8605 × 10 ⁻⁹	1.3136 × 10 ⁻⁷	1.2071 × 10 ⁻⁸	7.3910 × 10 ⁻⁸
Combined	0.9999999921	0.9999998686	0.999999879	0.9999999261



Troubleshooting

Observation	Possible cause	Recommended action
Faint or no signal from both the DNA Control 007 and the DNA	The incorrect volume of Master Mix or Primer Set was used.	Use the correct volume of Master Mix or Primer Set.
test samples at all loci, including the IQC markers	The DNA Polymerase was not activated.	Repeat the amplification with an initial hold at 95°C for 1 minute.
	The Master Mix was not vortexed thoroughly before aliquoting.	Vortex the Master Mix thoroughly.
	The Primer Set was exposed to too much light.	Replace the Primer Set and store it protected from light.
	Evaporation.	Ensure that the plate is properly sealed with film and that a compression pad was used with the GeneAmp™ PCR System 9700. (A compression pad should not be used with other validated thermal cyclers.)
	The thermal cycler malfunctioned.	See the thermal cycler user manual and check the instrument calibration.
	Incorrect thermal cycler conditions were used.	Use correct thermal cycler conditions.
	A MicroAmp [™] base was used with a tray/retainer set and tubes in GeneAmp [™] PCR System 9700.	Remove the MicroAmp [™] base.
	The tubes or plate were not seated tightly in the thermal cycler during amplification.	Push the tubes or plate firmly into the block after first cycle.
	The wrong PCR reaction tubes or plate were used.	Use MicroAmp [™] Reaction Tubes with Caps or the MicroAmp [™] Optical 96-well Reaction Plate for the GeneAmp [™] PCR System 9700 or Veriti [™] Thermal Cycler.
	Insufficient PCR product was electrokinetically injected.	Use correct genetic analyzer settings.
	Degraded formamide was used.	Check the storage of formamide. Do not thaw and refreeze multiple times. Try Hi-Di [™] Formamide.
Positive signal from DNA Control 007 but partial or no signal from DNA test samples when IQC peaks are present and balanced	The quantity of test DNA sample is below the assay sensitivity.	Quantify DNA and (when possible) add 500 pg of DNA. For low concentration samples, add up to 15 µL of the DNA sample to the reaction mix.

Observation	Possible cause	Recommended action
Positive signal from DNA Control 007 but partial or no signal from DNA test samples when IQC peaks are present and balanced	The test sample DNA is severely degraded.	Use the Quantifiler [™] HP DNA Quantification Kit or the Quantifiler [™] Trio DNA Quantification Kit to evaluate sample quality during the quantification step. If DNA is degraded, reamplify with an increased amount of DNA or consider using the Precision ID GlobalFiler [™] NGS STR Panel.
	The test sample was diluted in the wrong buffer (for example, a TE buffer with an incorrect EDTA concentration).	Redilute DNA using low-TE buffer (with 0.1 mM EDTA).
Positive signal from DNA Control 007 but partial or no signal from DNA test samples when IQC peaks are present	The test sample contains a high concentration of PCR inhibitor (for example, heme compounds, certain dyes).	Quantify the DNA, then use the minimum necessary volume of test sample DNA. Wash the sample in a Centricon [™] -100
and unbalanced	The test sample was diluted in the wrong buffer (for example, a TE buffer with an incorrect EDTA concentration).	centrifugal filter unit. Redilute DNA using low-TE buffer (with 0.1 mM EDTA).
Positive signal from DNA Control 007 and elevated signal from DNA test samples when IQC peaks are present and unbalanced	The quantity of the test sample DNA is in excess of recommended input amount, which can cause loss of balance in IQC peaks.	Quantify DNA, then use 500 pg.
More than two alleles present at a locus	Exogenous DNA is present in the sample.	Use appropriate techniques to avoid introducing foreign DNA during laboratory handling.
	Stutter product (–1 repeat unit position) was amplified.	Ensure that stutter filters are applied.
	Stutter filters apply to sample input amounts equal or greater than 250 pg. Samples well below the recommended input amount may exhibit stutters values exceeding the filters due to the stochastic effects of the PCR.	Increase the sample input above 250 pg, if possible. (Note, the optimum sample input is 500 pg.)
	The test sample contained mixed DNA.	If a mixed profile is not expected, check that laboratory protocols relating to cleanliness are followed.
	Incomplete 3´A base addition (n-1 nt position) occured.	Include the final extension step of 60°C for 5 minutes in the PCR.
		Remove amplified plate from storage (thaw if necessary) and place on thermal cycler at 60°C for 15 minutes.
		Check the quantity of the original sample DNA to ensure input is less than 750 pg per reaction. Adjust input as necessary during reamplification.

Appendix A Troubleshooting Population data

Observation	Possible cause	Recommended action
More than two alleles present at a locus		If the total amount of DNA in the reaction exceeds 1 ng, adjust the final extension time to 15 minutes to minimize incomplete 3' A base addition.
	The signal exceeds the dynamic range of the instrument and is causing signal "pull-up" into adjacent channels.	Check that you are using the recommended number of PCR cycles. Repeat PCR amplification using reduced input DNA amount, or interpret the off-scale data according to your laboratory procedure.
		Check that you are using the recommended injection conditions on the instrument.
	Poor spectral separation occurred.	Perform a spectral calibration.
		Confirm that Filter Set J6-T modules are installed and used for analysis.
	Too much DNA was present in the reaction.	Use the recommended amount of template DNA: 500 pg for 30 PCR cycles.
	The double-stranded DNA was not completely denatured.	Use the recommended amount of Hi-Di [™] Formamide and heat the sample plate at 95°C for 3 minutes.
Poor peak height balance	Incorrect thermal cycler conditions were used.	Use correct thermal cycler conditions.
Some but not all loci visible on electropherogram of DNA Test Samples	The PCR reaction volume you used is lower than the volume required for the amplification.	Use the correct PCR reaction volume: 25 µL
STR profiles contain many off- scale alleles	DNA quantification was not performed or not accurate.	Ensure that DNA quantification is accurate.



Materials required but not supplied

Unless otherwise indicated, all materials are available through **thermofisher.com**. MLS: Fisher Scientific (**fisherscientific.com**) or other major laboratory supplier.

STR kit required materials

Item	Source
NGM Detect [™] PCR Amplification Kit, 200 reactions	A31832

Sample preparation required materials

Item	Source
GeneScan [™] 600 LIZ [™] Size Standard v2.0, 2 × 200 μL	4408399
IMPORTANT! Do not use GeneScan [™] 350 ROX [™] , GeneScan [™] 500 ROX [™] , or GeneScan [™] 500 LIZ [™] Size Standards with this kit.	
Low-TE buffer (10 mM Tris, 0.1 mM EDTA, pH 8.0)	Teknova T0223
	or see "Prepare low-TE buffer" on page 19.
Hi-Di [™] Formamide, 25-mL	4311320

Thermal cycler required materials

$\mathsf{ProFlex}^{^{\mathsf{TM}}} \mathsf{PCR} \mathsf{System}$

Item	Source
ProFlex [™] 96-well PCR System	4484075
ProFlex [™] 2 × 96-well PCR System	4484076
ProFlex [™] 3 × 32-Well PCR System	4484073

$Veriti^{^{\mathrm{TM}}}$ Thermal Cycler

Item	Source
Veriti [™] 96-Well Thermal Cycler	4479071
(Optional) Tabletop centrifuge with 96-Well Plate Adapters	MLS

GeneAmp[™] PCR System 9700

Item	Source
GeneAmp [™] PCR System 9700, 96-Well Silver	N8050001
GeneAmp [™] PCR System 9700, 96-Well Gold-Plated	4314878
Silver 96-Well Sample Block	N8050251
Gold-Plated 96-Well Block	4314443

Genetic analyzer required materials

3500 Series Genetic Analyzer

Item	Source
3500 Series Data Collection Software 3.1 Upgrade (RUO)	A26287 ^[1]
3500 Series Data Collection Software 3.1 (RUO)	4475183 ^[1]
HID Updater 3500 Data Collection Software v2	4480670
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
DS-37 Matrix Standard Kit (Dye Set J6-T)	A31234
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

Item	Source
96-Well Septa for 3500/3500xL Genetic Analyzers	4412614
Septa Cathode Buffer Container, 3500 series	4410715

^[1] Contact your Thermo Fisher Scientific HID representative.

3130 Series Genetic Analyzer required materials

item	Source
3130 Data Collection Software v4	4475105
3130xl Data Collection Software-v4	4475126
3130/3730 Data Collection Software-v4 6-Dye Module v1	4478404
96-Well Plate Septa	4315933
Reservoir Septa	4315932
3130/3130 <i>xl</i> Genetic Analyzer 16-Capillary Array, 36 cm	4315931
POP-4 [™] Polymer for the 3130/3130 <i>xl</i> Genetic Analyzer	4352755
Running Buffer, 10×	402824
DS-37 Matrix Standard Kit (Dye Set J6-T)	A31234
MicroAmp [™] Optical 96-Well Reaction Plate	N8010560

Analysis software required materials

GeneMapper $^{\mathsf{TM}}$ *ID-X* Software v1.5.2 patch

In addition to the GeneMapper $^{\text{\tiny TM}}$ *ID-X* Software listed in the following table, a v.1.5.2 patch is required. The patch enables full functionality of the data quality assessment tools when using the NGM Detect $^{\text{\tiny TM}}$ PCR Amplification Kit.

The patch is available for free download at thermofisher.com/us/en/home/technical-resources/software-downloads/genemapper-id-x-software.html.

$GeneMapper^{TM}$ /D-X Software

Item	Source
GeneMapper [™] /D-XSoftware v1.5 Full Installation	A27884
GeneMapper [™] /D-XSoftware v1.5 Client Installation	A27886
GeneMapper [™] /D-XSoftware v1.4 Full Installation	4479707
GeneMapper [™] /D-XSoftware v1.4 Client Installation	4479711

Miscellaneous required materials

Plates and tubes

item	Source
MicroAmp [™] 96-Well Tray	N8010541
MicroAmp [™] Reaction Tube with Cap, 0.2 mL	N8010540
MicroAmp [™] 8-Tube Strip, 0.2 mL	N8010580
MicroAmp [™] Optical 8-Cap Strips	4323032
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

Laboratory supplies

Item	Source	
Various procedures		
Aerosol resistant pipette tips	MLS ^[1]	
Microcentrifuge tubes	MLS	
Pipettors	MLS	
Tape, labeling	MLS	
Tube, 50-mL Falcon [™]	MLS	
Tube decapper, autoclavable	MLS	
Deionized water, PCR grade	MLS	
Vortex	MLS	

^[1] Major laboratory supplier



PCR work areas

Work area setup and lab design	131
PCR setup work area materials	131
Amplified DNA work area	132

Work area setup and lab design

Many resources are available for the appropriate design of a PCR laboratory. If you are using this kit for:

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

The sensitivity of this 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).

Process samples carefully 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 materials

IMPORTANT! Do not remove these items from the PCR Setup Work Area.

- Calculator
- Gloves, disposable
- Marker pen, permanent
- Microcentrifuge
- Microcentrifuge tubes, 1.5-mL, or 2.0-mL, or other appropriate nuclease-free 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.

Use only the validated thermal cyclers listed in "Instruments and software compatibility" on page 15.



Safety



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, and so on). To obtain SDSs, see the "Documentation and Support" section in this document.

Appendix D Safety Chemical safety

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. 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! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological* and *Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:
 - www.cdc.gov/biosafety/publications/bmbl5/BMBL.pdf
- World Health Organization, Laboratory Biosafety Manual, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:
 - www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf

Documentation and support

Related documentation

Document title	Pub. No.
STR kits	
$NGM\ Detect^{^{ imes}}\ PCR\ Amplification\ Kit$ – $PCR\ Amplification\ and\ CE\ Quick\ Reference$	100044088
NGM Detect [™] PCR Amplification Kit – PCR Setup Quick Reference	100044087
Technical Note: Updated NGM Detect [™] PCR Amplification Kit: Validation and Comparative Study	_
Note: Go to thermofisher.com/document-connect/document-connect.html?url=https://assets.thermofisher.com/TFS-Assets/GSD/Technical-Notes/ngm-detect-results-tech-note.pdf to open the Technical Note.	
Quantification kits	
Quantifiler [™] HP and Quantifiler [™] Trio DNA Quantification Kits User Guide	4485355
Thermal cyclers	
ProFlex [™] PCR System User Guide	MAN0007697
Veriti [™] Thermal Cycler User Guide	4375799
GeneAmp [™] PCR System 9700 Base Module User Manual	4303481
3500 Series Genetic Analyzer	
3500/3500xL Genetic Analyzer with 3500 Series Data Collection Software v1 User Guide	4401661
3500/3500xL Genetic Analyzer with 3500 Series Data Collection Software v2 User Guide	4476988
HID Updater 3500 Data Collection Software v2.0 User Bulletin	NA
3500/3500xL Genetic Analyzer with 3500 Series Data Collection Software 3.1 User Guide	100031809
3500 Series Data Collection Software v3 User Bulletin: New Features and HID Validation Summary	MAN0010812
3500 Series Data Collection Software v3.1 User Bulletin: New Features and HID Validation Summary	MAN0014110
3130 <i>xl</i> Series Genetic Analyzer	
3130/3130xl Genetic Analyzers Maintenance, Troubleshooting, and Reference Guide	4352716
3130/3130xl Genetic Analyzers Using Data Collection Software v3.0 User Bulletin	4363787
3130/3130xl Genetic Analyzers Getting Started Guide	4352715

Document title	Pub. No.
3130/3130xl Genetic Analyzers Quick Reference Card	4362825
3130/3130xl Genetic Analyzers AB Navigator Software Administrator Guide	4359472
GeneMapper [™] <i>ID-X</i> Software all versions	
GeneMapper [™] ID-X Software Bin Overlap User Bulletin	100029546
GeneMapper [™] <i>ID-X</i> Software v1.0	
GeneMapper [™] ID-X Software v1.0 Getting Started Guide— Basic Features	4375574
GeneMapper [™] ID-X Software v1.0 Quick Reference— Basic Features	4375670
GeneMapper [™] ID-X Software v1.0 Installation Guide	4476603
GeneMapper [™] ID-X Software v1.0 Administrator Guide	4376327
GeneMapper [™] ID-X Software v1.0 Reference Guide	4375671
GeneMapper [™] <i>ID-X</i> Software v1.1	
GeneMapper [™] ID-X Software v1.1 Getting Started Guide— Mixture Analysis Tool	4396773
GeneMapper [™] <i>ID-X</i> Software v1.2	
GeneMapper [™] ID-X Software v1.2 Verification Experiments and Installation Procedures User Bulletin	4462639
GeneMapper [™] ID-X Software v1.2 Quick Reference— Mixture Analysis Tool	4426482
GeneMapper [™] ID-X Software v1.2 Reference Guide	4426481
GeneMapper [™] <i>ID-X</i> Software v1.3	
GeneMapper [™] ID-X Software v1.3 Verification Experiments and Installation Procedures User Bulletin	4470483
GeneMapper [™] / <i>D-X</i> Software v1.4	
GeneMapper [™] ID-X Software v1.4 New Features and Installation Procedures User Bulletin	4477684
GeneMapper [™] / <i>D-X</i> Software v1.5	
GeneMapper [™] ID-X Software v1.5 New Features and Verification User Bulletin	100031708
GeneMapper™ ID-X Software v1.5 Getting Started Guide— Basic Features	100031701
GeneMapper™ ID-X Software v1.5 Quick Reference— Basic Features	100031702
GeneMapper™ ID-X Software v1.5 Getting Started Guide— Mixture Analysis Tool	100031704
GeneMapper™ ID-X Software v1.5 Quick Reference— Mixture Analysis Tool	100031705
GeneMapper™ ID-X Software v1.5 Installation Guide	100031706
GeneMapper™ ID-X Software v1.5 Administrator Guide	100031703
GeneMapper [™] ID-X Software v1.5 Reference Guide	100031707

Customer and technical support

For support:

- In North America—Send an email to HIDTechSupport@thermofisher.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 **thermofisher.com/support** to obtain the following information.

- Worldwide contact telephone numbers
- Product support
- Order and web support
- Safety Data Sheets (SDSs; also known as MSDSs)

Additional product documentation, including user guides and Certificates of Analysis, are available by contacting Customer Support.

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 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.

References

Akane, A., Matsubara, K., Nakamura, H., Takahashi, S., and Kimura, K. 1994. Identification of the heme compound copurified with deoxyribonucleic acid (DNA) from bloodstains, a major inhibitor of polymerase chain reaction (PCR) amplification. *J. Forensic Sci.* 39:362–372.

Barber, M.D. 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., 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., 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.

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.* 148:249–258.

Bender, K., Farfan, M.J., Schneider, P.M. 2004. Preparation of degraded human DNA under controlled conditions. *Forensic Sci. Int.* 139:134–140.

Brinkmann, 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.

Budowle, B. et al. 2010. Population genetic analyses of the NGM STR loci. *Int. J. Legal Med.* e-publication **www.springerlink.com/content/p713q3w5440674u3/**

Butler, J.M. 2005. Forensic DNA Typing. Burlington, MA:Elsevier Academic Press.

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.

Clark J.M. 1988. Novel non-templated nucleotide addition reactions catalyzed by procaryotic and eucaryotic DNA polymerases. *Nucleic Acids Res.* 16:9677–9686.

DeFranchis, R., Cross, N.C.P., Foulkes, N.S., and Cox, T.M. 1988. A potent inhibitor of Taq DNA polymerase copurifies with human genomic DNA. *Nucleic Acids Res.* 16:10355. DNA Advisory Board, Federal Bureau of Investigation, U.S. Department of Justice. 1998. Quality assurance standards for forensic DNA testing laboratories.

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.

Excoffier, L., Lischer, H.E.L. 2010. A new series of programs to perform population genetics analyses under Linux and Windows. Arleguin suite v. 3.5. *Mol. Ecol. Res.* 10:564–567.

Frank, W., Llewellyn, B., Fish, P., et al. 2001. Validation of the AmpFlSTR Profiler Plus PCR Amplification Kit for use in forensic casework. *J. Forensic Sci.* 46:642–646.

Holt, C., Stauffer, C., Wallin, J., et al. 2000. Practical applications of genotypic Surveys for forensic STR testing. *Forensic Sci. Int.* 112:91–109.

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., Matise, T.C. 2004. A combined linkage-physical map of the human genome. *Am. J. Hum. Genet.* 75:1143–1148.

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.

Li, B., Ge, J., Wu, F., Ye, L., Budowle, B., Vhen, Y. 2013. Population genetic analyses of the STR loci of the AmpFlSTR NGM SElect PCR Amplification Kit for Han population in Fujian Province, China. *Int J. Legal Med.* 127:345–346.

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 nontemplated 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.

Möller, 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.

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.

Puers C., Hammond H.A., Jin L., Caskey C.T., Schumm J.W. 1993. 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. *Am J. Hum. Genet.* 53(4):953–958.

Raymond M. and Rousset F. 1995. GENEPOP (version 1.2): population genetics software for exact tests and ecumenicism. *J. Heredity* 86:248–249.

Rousset, F. 2008. Genepop'007: A complete reimplementation 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.

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.

Scientific Working Group on DNA Analysis Methods (SWGDAM). 2016. Validation Guidelines for DNA Analysis Methods. Available at http://media.wix.com/ugd/4344b0_813b241e8944497e99b9c45b163b76bd.pdf Accessed 22 February 2017.

Wallin, J.M., Buoncristiani, M.R., Lazaruk, K.D., Fildes, N., Holt, C.L., Walsh, P.S. 1998. SWGDAM validation of the AmpFlSTR 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.

Weir, B. 1990. Genetic Data Analysis. Sinauer Associates Sunderland, MA.

Index

+A 85 +A nucleotide addition 85 3' A 85, 86 3130 instrument 28 3500 instrument 24 6-dye license activation 28	references 24, 28 run module 24, 28 setup of the 3130 instrument 28 setup of the 3130xl instrument 28 setup of the 3500 and 3500xL instruments 24 extra peaks 77
spectral calibration 26, 30	GeneScan 600 LIZ Size Standard v2.0 48
600 LIZ Size Standard v2.0 48	GeneScan size standard, about 11
A	1
accuracy and reproducibility 63 alleles, off-ladder 65 allelic ladder, requirements for electrophoresis 23 artifacts 87	import panels, bins, and marker stutter 36 instrument and software compatibility 15
	L
В	limited product warranty 138 LIZ size standard
bins, import 36	about 11
biohazard safety 135	peak sizes 48 volume per reaction 27, 30 LIZ Size Standard v2.0 48
C	
characterization of loci, validation 88 control DNA	М
007 11	marker stutter, import 36
profile 13	materials not supplied 127
D	Р
developmental validation 57 direct amplification 22 DNA control profile 13 documentation, related 136 DS-37 matrix standard 26, 30 dye set for 6-dye samples 26, 30	panels check version 35 import 36 PCR conditions 21 perform 21 setup 131
E	work areas 131
electrophoresis data collection software 24, 28	Q
prepare samples 27, 30	quantification, DNA 17

R

required materials 127 run module for electrophoresis 3130 instrument 28 3130xl instrument 28 3500 and 3500xL instruments 24

S

safety, biohazard 135 sensitivity 92 size standard 48 spectral calibration 26, 30 stutter, peaks 83 stutter file, import 36

T

terms and conditions 138 thermal cyclers for use with kit 15 programming 21 troubleshooting 124

٧

validation, importance 56

W

warranty 138 work area, PCR setup 131

