

# Verifiler™ Plus PCR Amplification Kit

## USER GUIDE

**Catalog Number** A35495

**Publication Number** MAN0017493

**Revision** D

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

**ThermoFisher**  
S C I E N T I F I C



Revision history: MAN0017493 D (English)

Revision	Date	Description
D	22 April 2025	The following changes were made: <ul style="list-style-type: none"><li>• Storage conditions were updated (see “Contents and storage” on page 14).</li><li>• Compatible instruments, compatible software, and materials required were updated (throughout the document).</li><li>• Copy edits and formatting changes were made to align with current documentation style (throughout the document).</li></ul>
C.0	10 August 2020	Validation experiments were added for the updated formulation. Information about the D10S2148 marker was added.
B.0	17 October 2018	A new chapter was added: Chapter 6, “Experiments and results”. Minor edits.
A.0	24 May 2018	New document for the Verifiler™ Plus PCR Amplification Kit.

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.

**Important Licensing Information:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, 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.

©2018-2025 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

■ CHAPTER 1	Product information .....	7
	Product description .....	7
	Internal Quality Control System for PCR .....	7
	About the primers .....	8
	Dyes used in the kit .....	9
	Loci amplified by the kit .....	10
	Standards and controls that are required .....	11
	Contents and storage .....	14
	Required materials not supplied .....	15
	Instruments and software compatibility .....	15
	Workflow .....	17
■ CHAPTER 2	Perform PCR .....	18
	Validated DNA amounts and PCR cycles .....	18
	DNA quantification .....	18
	Importance of DNA quantification before STR analysis .....	18
	Effect of DNA quantity on results .....	19
	Methods of quantifying DNA .....	19
	Before you begin .....	20
	<i>(Optional)</i> Prepare low-TE buffer .....	20
	<i>(Before first use of the kit)</i> Thaw reagents .....	20
	Prepare the amplification kit reactions .....	21
	Perform PCR .....	22
	Direct amplification .....	23
■ CHAPTER 3	Perform electrophoresis .....	24
	Allelic ladder requirements for electrophoresis .....	24
	<i>(Before first use of the kit)</i> Set up the capillary electrophoresis instrument .....	25
	Data collection software setup .....	25
	Perform spectral calibration .....	26
	Prepare samples for electrophoresis and start the run .....	28

■	<b>CHAPTER 4</b>	<b>Analyze data with GeneMapper™ <i>ID-X</i> Software</b>	<b>29</b>
		Overview of GeneMapper™ <i>ID-X</i> Software	29
		Allelic ladder requirements for data analysis	30
		File names and versions used in this section	30
		Set up the GeneMapper™ <i>ID-X</i> Software for analysis (before first use of the kit)	31
		Workflow	31
		Check panel, bin, and stutter file versions on your computer	31
		<i>(If needed)</i> Download newer versions of panel, bin, and stutter files	32
		<i>(If needed)</i> Import panels, bins, and marker stutter	32
		<i>(Optional)</i> Define custom table or plot settings	36
		Create an analysis method	37
		Create an analysis method	37
		Enter Analysis Method settings	37
		<i>(If needed)</i> Create a size standard definition file	45
		About the GS600_LIZ_(60–460) size standard definition file	46
		Create a size standard definition file	46
		Analyze and edit sample files with GeneMapper™ <i>ID-X</i> Software	48
		Examine or edit a project	48
		For more information on using the GeneMapper™ <i>ID-X</i> Software	48
■	<b>CHAPTER 5</b>	<b>Assess the PCR reaction with the Internal Quality Control System</b>	<b>49</b>
		Overview of the Internal Quality Control system	49
		Evaluate the PCR reaction	49
		Balanced profile	51
		Ski slope profile with decreased IQCL peak height	51
		Ski slope profile with IQC peaks	52
		No sample peaks with IQC peaks	52
■	<b>CHAPTER 6</b>	<b>Experiments and results</b>	<b>53</b>
		Importance of validation	53
		Experiment conditions	53
		Repeated validation experiments using the updated primer set formulation	54
		Laboratory requirements for internal validation	54
		Developmental validation	54
		SWGDM guideline 2.2.1	54
		SWGDM guideline 3.9.2	54
		SWGDM guideline 3.9.6	55
		PCR components	55
		Thermal cycling temperatures	55
		PCR cycle number	56

Accuracy, precision, and reproducibility .....	57
SWGDM guideline 3.5 .....	57
Accuracy observation .....	57
Precision and size window description .....	60
Precision and size window observation .....	60
Extra peaks in the electropherogram .....	76
Causes of extra peaks .....	76
Extra peaks: Stutter .....	76
Extra peaks: Addition of 3' A nucleotide .....	82
Extra peaks: Artifacts .....	83
Characterization of loci .....	85
SWGDM guideline 3.1 .....	85
Loci in this kit .....	85
Nature of polymorphisms .....	85
Inheritance .....	86
Mapping .....	86
Genetic linkage .....	86
Species specificity .....	87
SWGDM Guideline 3.2 .....	87
Nonhuman study observation .....	87
Sensitivity .....	89
SWGDM guideline 3.3 .....	89
Sensitivity observation .....	89
Stability .....	90
SWGDM guideline 3.4 .....	90
Lack of amplification of some loci .....	90
Degraded DNA .....	91
Effect of inhibitors .....	91
Mixture studies .....	94
SWGDM guideline 3.8 .....	94
Mixture study overview .....	94
Mixture study observation .....	95
Resolution of genotypes in mixed samples .....	95
Limit of detection of the minor component .....	96
Population data .....	98
SWGDM guideline 3.7 .....	98
Loci in the kit .....	98
Population distribution .....	99
Population samples used in these studies .....	99
Concordance studies .....	100
Probability of identity .....	102
Probability of paternity exclusion .....	119

- **APPENDIX A** Troubleshooting ..... 120
- **APPENDIX B** Materials required but not supplied ..... 123
  - Sample preparation (PCR or CE) required materials ..... 123
  - Thermal cycler required materials ..... 123
    - ProFlex™ PCR System ..... 123
    - VeritiPro™ Thermal Cycler ..... 124
  - Capillary electrophoresis instrument required materials ..... 124
    - SeqStudio™ Flex Series Genetic Analyzer for Human Identification ..... 124
    - SeqStudio™ Genetic Analyzer for HID ..... 125
    - 3500 Series Genetic Analyzer for Human Identification ..... 125
  - Analysis software required materials ..... 126
    - GeneMapper™ *ID-X* Software ..... 126
  - Miscellaneous required materials ..... 126
    - Plates and tubes ..... 126
    - Laboratory supplies ..... 127
- **APPENDIX C** PCR work areas ..... 128
  - Work area setup and lab design ..... 128
  - PCR setup work area materials ..... 128
  - Amplified DNA work area ..... 129
- **APPENDIX D** Safety ..... 130
  - Chemical safety ..... 131
  - Biological hazard safety ..... 132
- **APPENDIX E** Documentation and support ..... 133
  - Related documentation ..... 133
  - Customer and technical support ..... 135
  - Limited product warranty ..... 135
- References ..... 136



# Product information

■ Product description .....	7
■ Contents and storage .....	14
■ Required materials not supplied .....	15
■ Instruments and software compatibility .....	15
■ Workflow .....	17

---

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

---

## Product description

The Verifiler™ Plus PCR Amplification Kit is a 6-dye, short tandem repeat (STR) multiplex assay for the amplification of human genomic DNA.

The kit amplifies the following:

- 23 autosomal STR loci: D3S1358, vWA, D16S539, CSF1PO, D6S1043, D8S1179, D21S11, D18S51, D5S818, D2S441, D19S433, FGA, D10S1248, D22S1045, D1S1656, D13S317, D7S820, Penta E, Penta D, TH01, D12S391, D2S1338, and TPOX
- 1 insertion/deletion polymorphic marker on the Y chromosome: Y indel
- Amelogenin (sex determining marker)
- 2 Internal Quality Control (IQC) markers; see “Internal Quality Control System for PCR” on page 7

## Internal Quality Control System for PCR

The Verifiler™ Plus kit includes an Internal Quality Control (IQC) System for PCR. The IQC System has two synthetic targets that are amplified with the sample:

- IQCS (small), a low molecular weight amplicon
- IQCL (large), a high molecular weight amplicon

The behavior of the IQC target peaks can be used to evaluate the success of the PCR reaction and to indicate sample quality. For more information, see Chapter 5, “Assess the PCR reaction with the Internal Quality Control System”.

## About the primers

The Verifiler™ Plus PCR Amplification Kit primers are manufactured to maximize the assay signal-to-noise ratio and simplify the interpretation of results.

For a comparison of the Verifiler™ Plus kit primers to other kit primers, see “Primer comparison” on page 8.

---

**Note:** The Verifiler™ Plus kit primers do not contain any mobility modifiers.

---

### Primer comparison

To compare the Verifiler™ Plus kit primers to the primers in the Verifiler™ Express PCR Amplification Kit, GlobalFiler™ PCR Amplification Kit, GlobalFiler™ Express PCR Amplification Kit, and NGM Detect™ PCR Amplification Kit, use Table 1. Do not use the table to make comparisons between the other kits; for example, do not compare the Verifiler™ Express kit to the GlobalFiler™ kit.

---

**Note:** A concordance population study was performed and genotypes were compared between the same samples amplified with the Verifiler™ Plus kit and the Verifiler™ Express kit and Huaxia™ Platinum™ kit. For more information, see “Concordance studies” on page 100.

---

**Table 1** Verifiler™ Plus kit primer comparison

Marker	Verifiler™ Plus kit comparison		
	Verifiler™ Express kit and Huaxia™ Platinum™ kit	GlobalFiler™ kit and GlobalFiler™ Express kit	NGM Detect™ kit
IQCS	Marker not included	Marker not included	Primers are the same
IQCL	Marker not included	Marker not included	Primers are the same
D2S441	1 primer is the same	1 primer is the same	Primers are different
D10S1248	Primers are different	Primers are different	Primers are different
D22S1045 <sup>[1]</sup>	Primers are the same	Primers are the same	Primers are different
D3S1358 <sup>[1]</sup>	Primers are different	Primers are different	Primers are different
D8S1179 <sup>[1]</sup>	Primers are the same	Primers are the same	Primers are different
D19S433 <sup>[1]</sup>	1 primer is the same	1 primer is the same	Primers are different
D5S818	Primers are different	Primers are different	Marker not included
vWA <sup>[1]</sup>	Primers are the same	Primers are the same	Primers are different
D1S1656	Primers are different	Primers are different	Primers are different
TH01	1 primer is the same	1 primer is the same	Primers are different
D21S11	Primers are the same	Primers are the same	Primers are different
D13S317	1 primer is the same	1 primer is the same	Marker not included



Table 1 Verifiler Plus kit primer comparison (continued)

Marker	Verifiler™ Plus kit comparison		
	Verifiler™ Express kit and Huaxia™ Platinum™ kit	GlobalFiler™ kit and GlobalFiler™ Express kit	NGM Detect™ kit
D12S391	Primers are different	Primers are different	Primers are different
D16S539 <sup>[1]</sup>	1 primer is different	1 primer is different	Primers are different
FGA <sup>[1]</sup>	Primers are the same	Primers are the same	Primers are different
D18S51	1 primer is the same	1 primer is the same	Primers are different
D7S820	Primers are different	Primers are different	Marker not included
D2S1338	Primers are the same	Primers are the same	Primers are different
CSF1PO	1 primer is the same	1 primer is the same	Marker not included
TPOX	1 primer is the same	1 primer is the same	Marker not included
D6S1043 <sup>[1]</sup>	Primers are different	Marker not included	Marker not included
Penta D <sup>[1]</sup>	Primers are different	Marker not included	Marker not included
Penta E	1 primer is the same	Marker not included	Marker not included
Y indel	Primers are the same	Primers are the same	Primers are the same
AMEL <sup>[1]</sup>	Primers are the same	Primers are the same	Primers are the same

<sup>[1]</sup> The Verifiler™ Plus kit has one or more additional SNP-specific primers.

## 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

## Loci amplified by the kit

Locus designation	Chromosome location	Alleles in the allelic ladder	Dye label	Alleles in DNA Control 007
IQCS	—	1, 2	6-FAM™	2
D3S1358	3p21.31	9–20		15, 16
vWA	12p13.31	11–24		14, 16
D16S539	16q24.1	5, 8–15		9, 10
CSF1PO	5q33.3–34	6–15		11, 12
D6S1043	6q15	7–25		12, 14
IQCL	—	1, 2		2
Y indel	Yq11.221	1, 2	VIC™	2
Amelogenin	X p22.1–22.3, Y: p11.2	X, Y		X, Y
D8S1179	8q24.13	5–19		12, 13
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
D18S51	18q21.33	7, 9, 10, 10.2, 11–13, 13.2, 14, 14.2, 15–27		12, 15
D5S818	5q21–31	7–18		11
D2S441	2p14	8–11, 11.3, 12–17	TED™	14, 15
D19S433	19q12	5–12, 12.2, 13, 13.2, 14, 14.2, 15, 15.2, 16, 16.2, 17, 17.2, 18.2, 19.2		14, 15
FGA	4q28	13–26, 26.2, 27–30, 30.2, 31.2, 32.2, 33.2, 42.2, 43.2, 44.2, 45.2, 46.2, 47.2, 48.2, 49.2, 50.2, 51.2		24, 26
D10S1248	10q26.3	8–19		12, 15
D22S1045	22q12.3	8–19	TAZ™	11, 16
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
D13S317	13q22–31	5–16		11
D7S820	7q11.21–22	6–15		7, 12
Penta E	15q26.2	5–26		7, 12
Penta D	21q22.3	2.2, 3.2, 5–17	SID™	11, 12

(continued)

Locus designation	Chromosome location	Alleles in the allelic ladder	Dye label	Alleles in DNA Control 007
TH01	11p15.5	4–9, 9.3, 10–12, 13.3	SID™	7, 9.3
D12S391	12p13.2	14–19, 19.3, 20–27		18, 19
D2S1338	2q35–37.1	11–28		20, 23
TPOX	2p23–2per	5–15		8

## Standards and controls that are required

The Verifiler™ Plus kit requires the following standards and controls for PCR amplification, PCR product sizing, and genotyping:

Item	Description	Included in the kit
DNA Control 007	Positive control. Used to evaluate amplification efficiency and to evaluate STR genotyping using the kit allelic ladder. See “DNA Control 007 profile” on page 12.	Yes
Verifiler™ Plus Allelic Ladder	Developed for accurate characterization of the alleles amplified in the kit. The allelic ladder allows automatic genotyping of most of the reported alleles for the loci in the kit. See “Loci amplified by the kit” on page 10 and Figure 2.	Yes
GeneScan™ 600 LIZ™ Size Standard v2.0 (Cat. No. 4408399)	Used for obtaining sizing results. This standard, which has been evaluated as an internal size standard, yields precise sizing results for PCR products.	No (order separately)

## DNA Control 007 profile

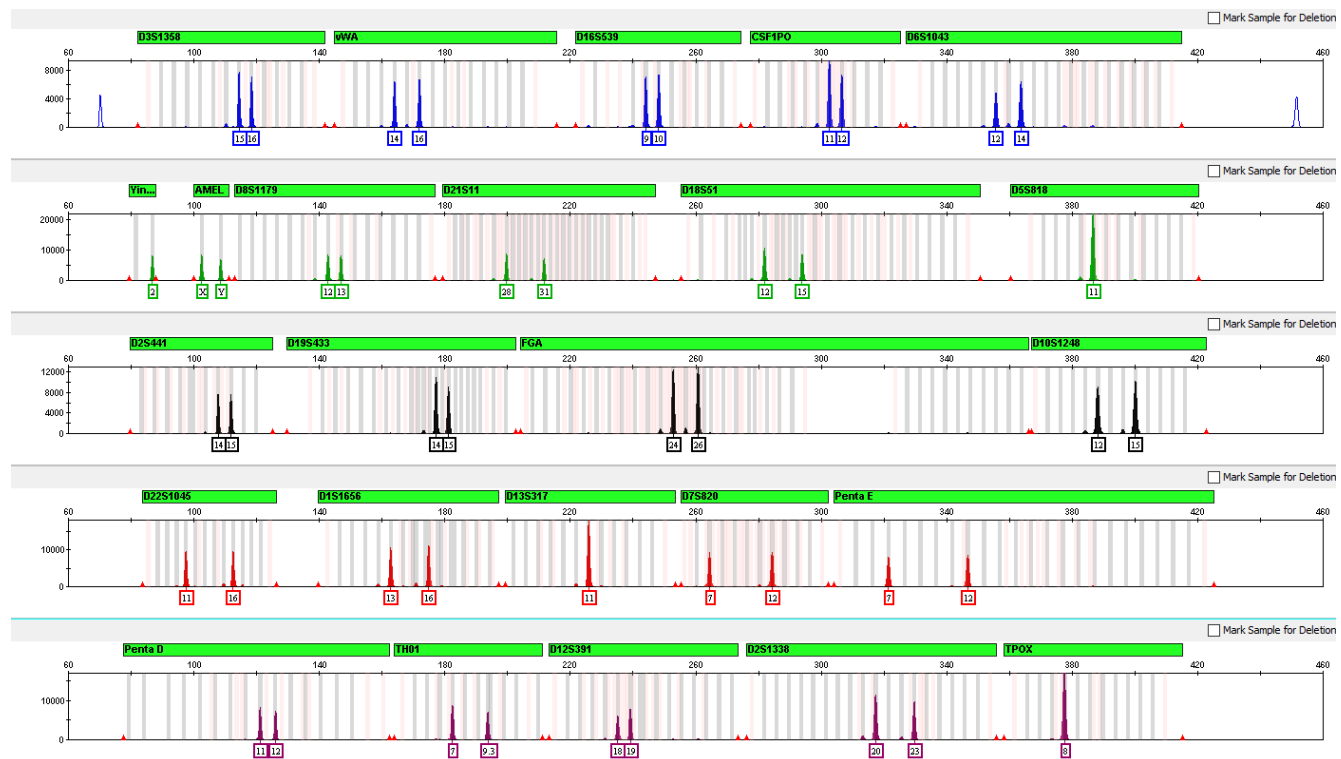
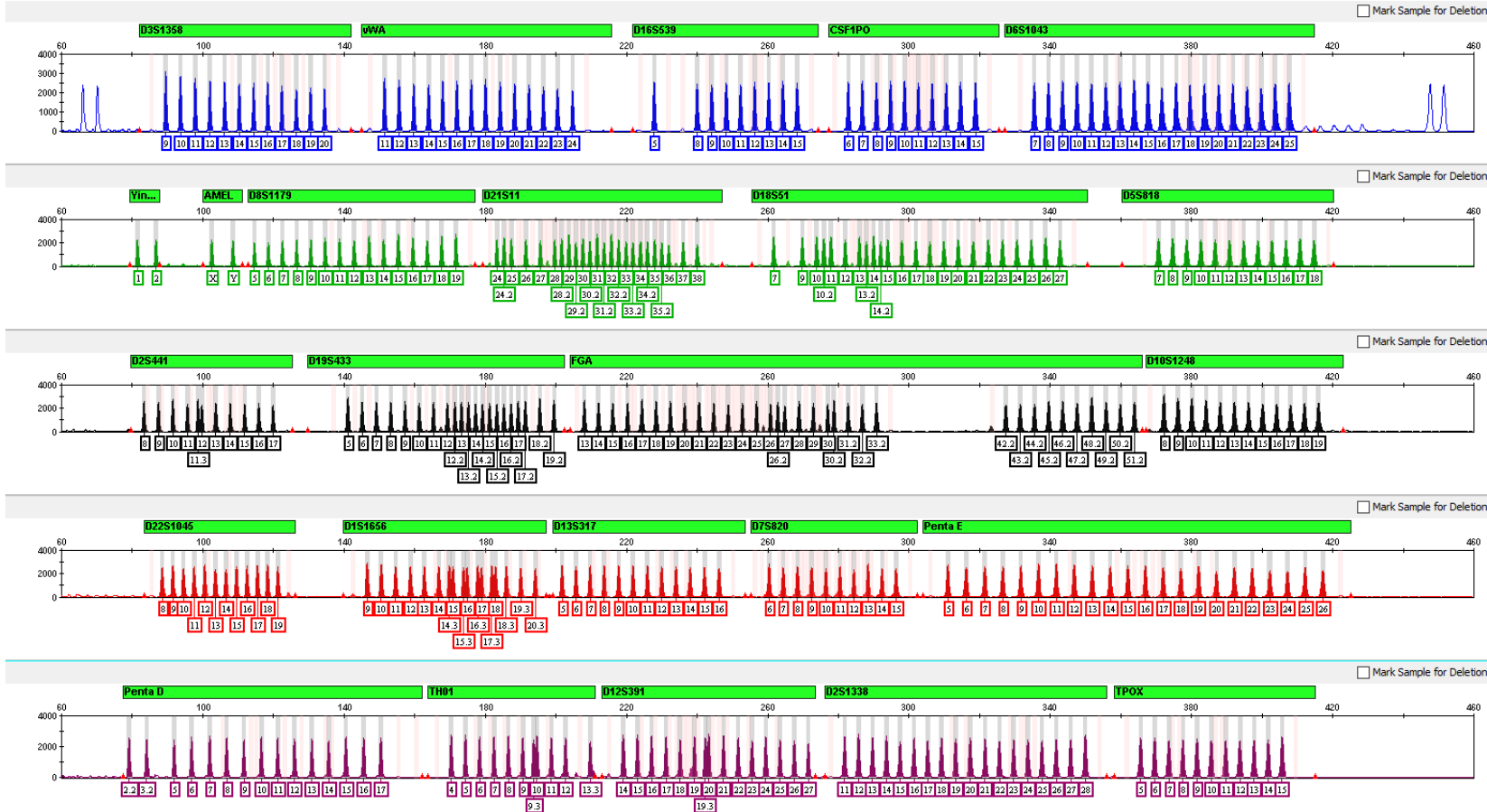


Figure 1 DNA Control 007 (500 pg) amplified with the Verifiler™ Plus kit and analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0–18,000 RFU) (updated formulation) The smallest and largest peaks in the FAM™ dye (blue) channel are the IQCS and IQCL markers, respectively.

## Allelic ladder profile



**Figure 2 GeneMapper™ ID-X Software plot of the Verifiler™ Plus Allelic Ladder (updated formulation)** The IQCS and IQCL markers are visible as the smallest and largest markers in the FAM™ dye (blue) channel. Each IQC marker has two alleles in the allelic ladder (alleles 1 and 2). However, the IQC markers are not normally analyzed and genotyped because they are primarily intended to be qualitative indicators.



## Contents and storage

The Verifiler™ Plus kit (Cat. No. [A35495](#)) contains sufficient quantities of the reagents for 200 amplification reactions at 25 µL/reaction.

---

**IMPORTANT!** The fluorescent dyes attached to the primers are light-sensitive. Protect the primer set and allelic ladder from light when not in use.

---



---

**IMPORTANT!** The allelic ladder contains PCR products that should not be amplified. To avoid contamination, store the allelic ladder separately from the other kit components and unamplified DNA.

---



---

**Note:** If there is more than one tube or bottle for a single reagent, thaw only the number of tubes or bottles required for the current number of reactions.

---



---

**IMPORTANT!** Do not refreeze kit components after thawing.

---



---

**Note:** For recommendations on kit handling, go to [thermofisher.com](http://thermofisher.com), then search for *Technical Note: Handling STR Kits and Ladder Decontamination*, or contact your local Human Identification representative.

---

Item	Description	Amount	Storage <sup>[1]</sup>
		200 reactions (Cat. No. <a href="#">A35495</a> )	
Verifiler™ Plus Master Mix	Contains enzyme, salts, dNTPs, bovine serum albumin, and 0.05% sodium azide in buffer and salt.	2 × 0.50 mL	–25°C to –15°C on receipt. 2–8°C after first use, up to the expiration date stated on the kit.
DNA Control 007	Contains 0.1 ng/µL human male genomic DNA from cell line in 0.05% sodium azide and buffer. <sup>[2]</sup> See “DNA Control 007 profile” on page 12 for information.	1 × 0.3 mL	
Verifiler™ Plus Primer Set	Contains forward and reverse primers to amplify human DNA targets.	2 × 0.25 mL	–25°C to –15°C on receipt. 2–8°C after first use, up to the expiration date stated on the kit.
Verifiler™ Plus Allelic Ladder	Contains amplified alleles. See Figure 2.	1 × 0.065 mL	Store protected from light.

<sup>[1]</sup> See packaging for expiration date. Do not use expired product.

<sup>[2]</sup> DNA Control 007 is included at a concentration that is appropriate for use as an amplification control (that is, to provide confirmation that the kit reagents can 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

**Note:** Compatible instruments and software that have been discontinued are not listed in this user guide.

### Thermal cyclers

- HID VeritiPro™ Thermal Cycler, 96-well
- ProFlex™ 96-well PCR System
- ProFlex™ 2 × 96-well PCR System
- ProFlex™ 3 × 32-Well PCR System

### Genetic analyzers and data collection software

Genetic analyzer	Data collection software
SeqStudio™ Flex Series Genetic Analyzer for Human Identification	SeqStudio™ Flex Series Instrument Software v1.1.1
SeqStudio™ Genetic Analyzer for HID	SeqStudio™ Data Collection Software v1.2.5
	SeqStudio™ Data Collection Software v1.2.4
	SeqStudio™ Data Collection Software v1.2.1
3500 Series Genetic Analyzer for Human Identification	3500 Series HID Data Collection Software v4.0.1
	3500 Series Data Collection Software 4 (Windows™ 10 operating system)

### Analysis software

Genetic analyzer	Analysis software
SeqStudio™ Flex Series Genetic Analyzer for Human Identification	GeneMapper™ ID-X Software v1.7.2 or later
SeqStudio™ Genetic Analyzer for HID	GeneMapper™ ID-X Software v1.6 or later
3500 Series Genetic Analyzer for Human Identification	GeneMapper™ ID-X Software v1.5 or later

### **For more information**

- For the instruments and software used during the kit validation, see Chapter 6, “Experiments and results”.
- For testing information on specific platforms, see the instrument or software user documentation.
- For ordering information, see Appendix B, “Materials required but not supplied”.



## Workflow

### Verifiler™ Plus PCR Amplification Kit

#### Extract and quantify DNA

1. Extract DNA—Go to: [www.thermofisher.com/hid-sampleprep](http://www.thermofisher.com/hid-sampleprep)
2. Quantify DNA—See “DNA quantification” on page 18

#### Perform PCR

1. “Prepare the amplification kit reactions” on page 21
2. “Perform PCR” on page 22

#### Perform capillary electrophoresis

1. “(Before first use of the kit) Set up the capillary electrophoresis instrument” on page 25
2. “Prepare samples for electrophoresis and start the run” on page 28)

#### Analyze data

1. “Set up the GeneMapper™ ID-X Software for analysis (before first use of the kit)” on page 31
2. “Create an analysis method” on page 37
3. “(If needed) Create a size standard definition file” on page 45
4. “Analyze and edit sample files with GeneMapper™ ID-X Software” on page 48
5. “Examine or edit a project” on page 48
6. Chapter 5, “Assess the PCR reaction with the Internal Quality Control System”

# 2

## Perform PCR

- Validated DNA amounts and PCR cycles ..... 18
- DNA quantification ..... 18
- Before you begin ..... 20
- Prepare the amplification kit reactions ..... 21
- Perform PCR ..... 22
- Direct amplification ..... 23

### Validated DNA amounts and PCR cycles

The kit is validated for use with 500 pg of DNA for 29 PCR cycles. The DNA input volume is 17.5 µL, for a total reaction volume of 25 µL.

### DNA quantification

#### Importance of DNA quantification before STR analysis

DNA quantification can be used to determine the following:

- If the sample contains sufficient human DNA and/or human male DNA to proceed with short tandem repeat (STR) amplification.
- (When using the Quantifiler™ Trio DNA Quantification Kit) The relative quantities of human male and female DNA in a sample. Relative quantities can help you select the appropriate STR chemistry.
- The amount of sample to use in STR analysis applications.
- If PCR inhibitors are present in a sample. If inhibitors are present, the sample may require additional purification before proceeding to STR analysis.
- The DNA quality, in regards to the inhibition level and the DNA degradation level. DNA quality can help you determine the likelihood of recovery of STR loci with larger amplicon sizes.

---

**Note:** Highly degraded samples that cannot be recovered by STR analysis with capillary electrophoresis can be analyzed with the Precision ID NGS System and Panels. Optimized for degraded samples, the Precision ID Identity Panel enables discrimination of individuals similar to STR genotype match probabilities. The Precision ID Ancestry Panel infers biogeographical ancestry for investigative leads.

---

## 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 the following:

- Fluorescence intensity that exceeds the linear dynamic range for detection by the capillary electrophoresis instrument (“off-scale” data). 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 is 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 reaction could be extremely low. Unbalanced amplification of the alleles can occur because of stochastic fluctuation.

## Methods of quantifying DNA

Kit	Detects	How it works
Quantifiler™ HP DNA Quantification Kit (Cat. No. 4482911)	<ul style="list-style-type: none"> <li>• Total human DNA (two targets—one small amplicon and one larger amplicon)</li> <li>• Degraded DNA</li> </ul>	<ul style="list-style-type: none"> <li>• Uses 5' nuclease assays with multiple-copy target loci, for improved detection sensitivity:<sup>[1]</sup> <ul style="list-style-type: none"> <li>– The human-specific target loci are multiple-copy, and dispersed on various autosomal chromosomes</li> <li>– The primary quantification targets have relatively short amplicons (75–80 bases), to improve the detection of degraded DNA samples</li> </ul> </li> </ul>
Quantifiler™ Trio DNA Quantification Kit (Cat. No. 4482910)	<ul style="list-style-type: none"> <li>• Total human DNA (two targets—one small amplicon and one larger amplicon)</li> <li>• Human male DNA</li> <li>• Degraded DNA</li> </ul>	<ul style="list-style-type: none"> <li>• Uses features that maximize consistency of quantification:           <ul style="list-style-type: none"> <li>– Genomic targets have conserved primer- and probe-binding sites</li> <li>– Minimal copy number variation between different individuals and population groups</li> </ul> </li> <li>• Contains a Large Autosomal target with a longer amplicon (&gt;200 bases) to help determine if a DNA sample is degraded</li> </ul>

<sup>[1]</sup> The detection sensitivity of the Quantifiler™ HP Kit and the Quantifiler™ Trio kit is improved over the Quantifiler™ Duo Kit.

---

**Note:** For information on the Quantifiler™ kits, see the *Quantifiler™ HP and Quantifiler™ Trio DNA Quantification Kits User Guide* (Pub. No. 4485354).

---

## Before you begin

### (Optional) Prepare low-TE buffer

We recommend Nuclease-Free Water (not DEPC-Treated) (Cat. No. [AM9937](#)) for sample preparation. However, you can also use low-TE buffer. Prepare the low-TE buffer as described in this procedure or use TE Buffer (Cat. No. [12090015](#)).

---

**IMPORTANT!** Ensure that the low-TE buffer has final concentrations of 10 mM Tris-HCl and 0.1 mM EDTA, pH 8.0. Do not use regular TE buffer because it has a higher EDTA concentration that can chelate Mg<sup>2+</sup> ions, therefore reducing PCR amplification efficiency and robustness.

---

1. Mix the buffer components together.
  - 10 mL of 1 M Tris-HCl, pH 8.0
  - 0.2 mL of 0.5 M EDTA, pH 8.0
  - 990 mL of nuclease-free water

---

**Note:** Adjust the volumes proportionally for specific needs.

---

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

### (Before first use of the kit) Thaw reagents

Thaw the master mix and primer set.

---

**IMPORTANT!** The fluorescent dyes attached to the primers are light-sensitive. Protect the primer set and allelic ladder from light when not in use.

---

---

**IMPORTANT!** Thawing is required only before first use of the kit. After first use, the reagents are stored at 2–8°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 and allelic ladder from light when not in use.

1. Vortex the master mix and primer set for 3 seconds. Before opening the tubes or bottles, remove droplets from the caps by briefly centrifuging the tubes or tapping the bottles on the bench.
2. Pipet the required component volumes into an appropriately sized, clear (non-colored), polypropylene tube.

Component	Amount per reaction
Master mix	5.0 $\mu\text{L}$
Primer set	2.5 $\mu\text{L}$

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

3. Vortex the reaction mix for 3 seconds, then briefly centrifuge.
4. Pipet 7.5  $\mu\text{L}$  of the reaction mix into each well of a MicroAmp™ Optical 96-Well Reaction Plate or each MicroAmp™ tube.
5. *(If needed)* Adjust the sample input amount and volume.
  - If the total sample input amount is >500 pg of DNA, dilute with nuclease-free water or low-TE buffer to obtain a 500-pg input in a 17.5- $\mu\text{L}$  volume.
  - If the total sample input volume is <17.5  $\mu\text{L}$ , bring to volume with nuclease-free water or low-TE buffer to obtain a 17.5- $\mu\text{L}$  input volume.
6. Prepare the samples and controls as shown in the following table, then add to the appropriate wells of a MicroAmp™ Optical 96-Well Reaction Plate or to each MicroAmp™ tube.

Component	Amount per reaction
Negative control	17.5 $\mu\text{L}$ of nuclease-free water or low-TE buffer
Test sample	17.5 $\mu\text{L}$ of DNA
Positive control	Combine, then add to the reaction well or tube: <ul style="list-style-type: none"> <li>• 5 <math>\mu\text{L}</math> of DNA Control 007 (0.1 ng/<math>\mu\text{L}</math>)</li> <li>• 12.5 <math>\mu\text{L}</math> of nuclease-free water or low-TE buffer</li> </ul>

The final reaction volume (sample or control plus reaction mix) is 25  $\mu\text{L}$ .

7. Seal the plate with MicroAmp™ Clear Adhesive Film or MicroAmp™ Optical Adhesive Film, or cap the tubes.

---

**IMPORTANT!** We recommend adhesive film for plate sealing to provide a consistent seal across all wells and prevent evaporation. Do not use caps for the plate, which may not provide a consistent seal across all wells.

---

8. Vortex the plate or tubes at medium speed for 3 seconds.
9. Centrifuge the tubes or plate at 3,000 × g for ~30 seconds in a tabletop centrifuge (with plate holders, if using 96-well plates).

Proceed to “Perform PCR” on page 22.

## Perform PCR

---

**Note:** This protocol is for casework only. For a direct amplification protocol, go to [thermofisher.com](http://thermofisher.com), then search for the technical note *Direct Amplification of Reference Samples Using the Verifiler™ Plus PCR Amplification Kit*, or contact your local Human Identification representative.

---

**IMPORTANT!** The kit is optimized for use with the thermal cyclers that are listed in “Instruments and software compatibility” on page 15.

---

1. Program the thermal cycler.
  - a. Set the ramping mode to **9700 Simulation**.
  - b. Set the thermal cycling conditions as shown in the following table.

Initial incubation step	First stage (2 cycles)		Second stage (27 cycles)		Final extension	Final hold
	Denature	Anneal/extend	Denature	Anneal/extend		
HOLD	CYCLE (29 cycles) <sup>[1]</sup>				HOLD	HOLD
95°C 1 minute	96°C 10 seconds	62°C 90 seconds	96°C 10 seconds	59°C 90 seconds	60°C 5 minutes	4°C ≤24 hours <sup>[2]</sup>

<sup>[1]</sup> See “Validated DNA amounts and PCR cycles” on page 18.

<sup>[2]</sup> The infinity (∞) setting allows an unlimited hold time.

2. Load the plate or tubes into the thermal cycler, close the heated cover, then start the run.
3. When the run is complete, store the amplified DNA.

Storage time	Temperature
<2 weeks	2–8°C
>2 weeks	–25°C to –15°C

---

**IMPORTANT!** Protect the amplified DNA from light.

---

## Direct amplification

The Verifiler™ Plus PCR Amplification Kit is designed and optimized for casework sample processing (including a pre-amplification extraction and cleanup step) with Verifiler™ Express PCR Amplification Kit as the partner kit. The Verifiler™ Express kit is designed and optimized to perform direct amplification of reference samples.

Some laboratories have expressed a preference for a single PCR amplification kit capable of processing both casework and reference samples. To support those laboratories, we have optimized a direct amplification PCR protocol. We performed studies to demonstrate that the Verifiler™ Plus kit can be used for direct amplification of reference samples. Specifically, we evaluated the performance of the kit using a direct amplification of blood and buccal samples collected on FTA™ cards, NUCLEIC-CARD™ devices, Bode Buccal DNA Collectors, blood stain cards made out of filter paper, and two types of swabs (Puritan™ Cotton Swabs and 4N6FLOQSwabs™).

---

**Note:** The direct amplification uses a different PCR protocol. Perform validation studies before using this kit for direct amplification of single-source samples, or use the Verifiler™ Express kit. At PCR cycle numbers <29, the IQC peaks may not be detected.

---

For more information on direct amplification studies, go to [thermofisher.com](https://www.thermofisher.com), then search for the *Technical Note: Direct Amplification of Reference Samples Using the Verifiler™ Plus PCR Amplification Kit*, or contact your local Human Identification representative.

# 3

## Perform electrophoresis

- Allelic ladder requirements for electrophoresis ..... 24
- (Before first use of the kit) Set up the capillary electrophoresis instrument ..... 25
- Prepare samples for electrophoresis and start the run ..... 28

### 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
SeqStudio™ 24 Flex Genetic Analyzer	1 per injection	24 samples	23 samples + 1 allelic ladder
SeqStudio™ 8 Flex Genetic Analyzer	1 per 3 injections	8 samples	23 samples + 1 allelic ladder
SeqStudio™ Genetic Analyzer	1 per 6 injections	4 samples	23 samples + 1 allelic ladder
3500xL Genetic Analyzer	1 per injection	24 samples	23 samples + 1 allelic ladder
3500 Genetic Analyzer	1 per 3 injections	8 samples	23 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. To facilitate accurate genotyping of all samples in your laboratory environment, perform internal validation studies to verify the required allelic ladder injection frequency.

---

It is critical to genotype using an allelic ladder that is 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.



## (Before first use of the kit) Set up the capillary electrophoresis instrument

### Data collection software setup

To analyze PCR products generated by the kit, you can use the data collection software and run parameters provided in this section. See the appropriate table for your instrument.

**Note:** With 0.5 ng of input DNA, our studies indicate that the injection conditions provided in this section produce well-balanced profiles 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 system. Large deviations from the recommended injection parameters can affect the performance of the size standard and allelic ladder, therefore validation is recommended.

**Note:** For detailed procedures, see the appropriate user documentation for your instrument.

**Table 2 Software setup: SeqStudio™ Flex Series Genetic Analyzer for Human Identification**

SeqStudio™ Flex Data Collection Software	(Optional) Additional software	Run parameters
v1.1.1	<ul style="list-style-type: none"> <li>SAE Administrator Console v2.1</li> <li>SeqStudio™ Plate Manager Software v2.1, v2.1.1</li> <li>SeqStudio™ Flex Remote Monitoring Software</li> </ul>	<b>Injection protocol:</b> HID_Protocol_J6T_36_POP4(xl)
		<b>Size standard:</b> GS600 LIZ (60–460)
		<b>Dye set:</b> J6-T (DS-37)
		<b>Run module:</b> HID_J6T_36_POP4(xl)
		<b>Injection conditions:</b> 1.2 kV/15 seconds (xl: 24 seconds)
		<b>Run conditions:</b> 13 kV/1,550 seconds

**Table 3 Software setup: SeqStudio™ Genetic Analyzer for HID**

SeqStudio™ Data Collection Software	(Optional) Additional software	Run parameters	Plate setup
v1.2.1, v1.2.4, v1.2.5	<ul style="list-style-type: none"> <li>SAE Administrator Console v2.0, v2.1</li> <li>SeqStudio™ Plate Manager Software v1.2, v1.3</li> </ul>	<b>Run Module:</b> HID Analysis	<b>Kit:</b> Verifiler™ Plus kit
		<b>Injection conditions:</b> 1.2 kV/10 seconds	<b>Dye set:</b> J6-T (DS-37)
		<b>Run conditions:</b> 11 kV/1,120 seconds	<b>Size standard:</b> GS600 LIZ (60–460)

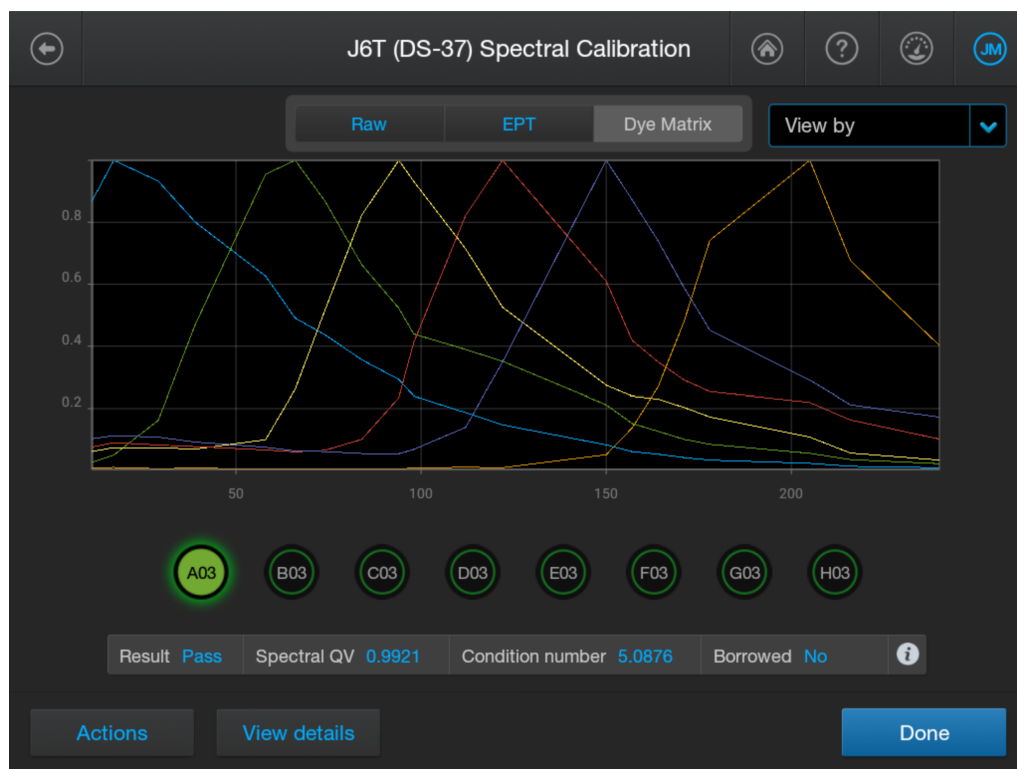
**Table 4 Software setup: 3500 Series Genetic Analyzer for Human Identification**

Operating system	3500 Data Collection Software	Run parameters
Windows™ 10	v4, v4.0.1	Assay: AB_J6T_LS_POP4(xl)
		Instrument protocol: AB_HID36_POP4(xl)_J6T_NT3200
		Run module: HID36_POP4(xl)
		Injection conditions: 1.2 kV/15 seconds (xl: 24 seconds)
		Run conditions: 13 kV/1,550 seconds
		Dye set: J6-T

## Perform spectral calibration

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

Examples of spectral calibrations are shown in this section. See the appropriate figure for your instrument.



**Figure 3 Example spectral calibration: SeqStudio™ Flex Series Genetic Analyzer for Human Identification**

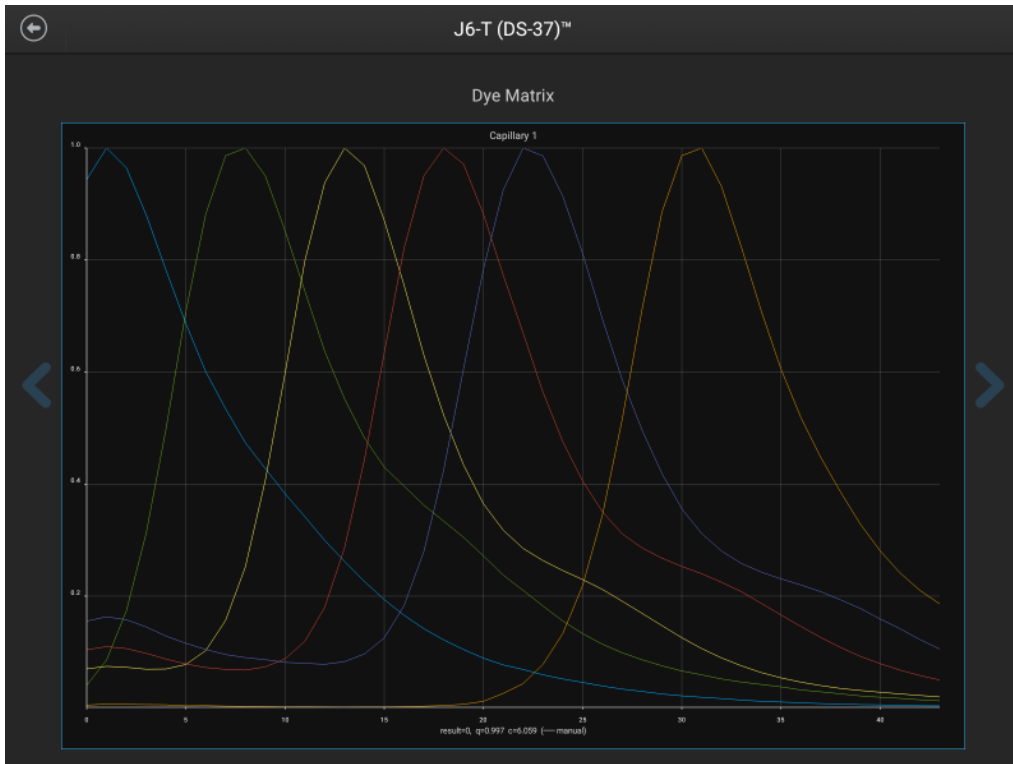


Figure 4 Factory-provided spectral calibration: SeqStudio™ Genetic Analyzer for HID

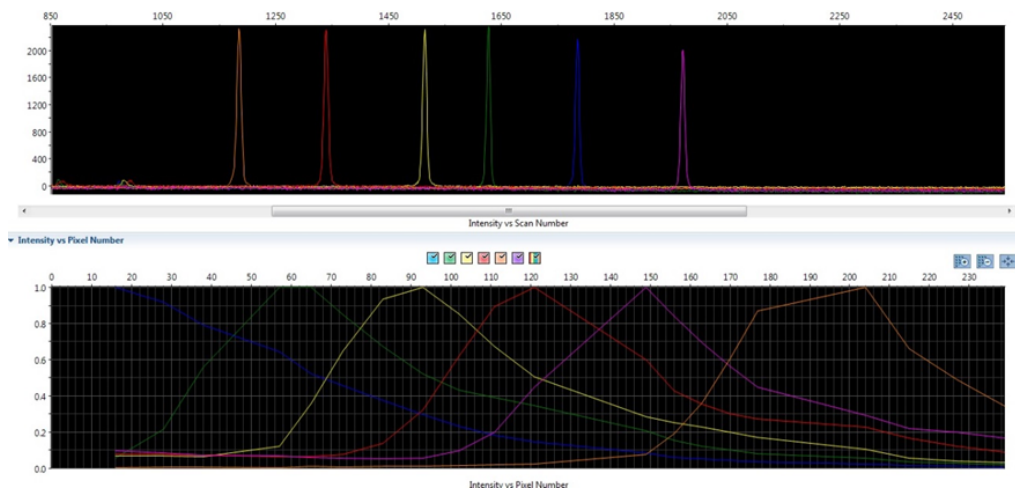


Figure 5 Example spectral calibration: 3500 Series Genetic Analyzer for Human Identification

## Prepare samples for electrophoresis and start the run

Prepare the samples for electrophoresis immediately before loading.

---

**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.

---

1. Pipet the required component amounts into an appropriately sized polypropylene tube.

Component	Amount per reaction
GeneScan™ 600 LIZ™ Size Standard v2.0	0.4 µL
Hi-Di™ Formamide	9.6 µL

---

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

---

**IMPORTANT!** The amount 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. Pipet the required component amounts into each well of a MicroAmp™ Optical 96-Well Reaction Plate.

Component	Amount per reaction
Formamide/size standard mixture	10 µL
PCR product or allelic ladder	1 µL

---

**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 for 3 minutes at 95°C.
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™ <i>ID-X</i> Software .....	29
■ Allelic ladder requirements for data analysis .....	30
■ File names and versions used in this section .....	30
■ Set up the GeneMapper™ <i>ID-X</i> Software for analysis (before first use of the kit) .....	31
■ Create an analysis method .....	37
■ (If needed) Create a size standard definition file .....	45
■ Analyze and edit sample files with GeneMapper™ <i>ID-X</i> Software .....	48
■ Examine or edit a project .....	48
■ For more information on using the GeneMapper™ <i>ID-X</i> Software .....	48

## Overview of GeneMapper™ *ID-X* Software

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

After capillary electrophoresis, the data collection software stores information for each sample in a FSA or HID file. The GeneMapper™ *ID-X* Software allows you to analyze and interpret the data from the FSA or HID files.

---

**Note:** For a list of GeneMapper™ *ID-X* Software versions that are compatible with your kit and capillary electrophoresis instrument, see “Instruments and software compatibility” on page 15.

---

## 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™ 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 help 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](https://www.thermofisher.com/support).

## Set up the GeneMapper™ ID-X Software for analysis (before first use of the kit)

### Workflow

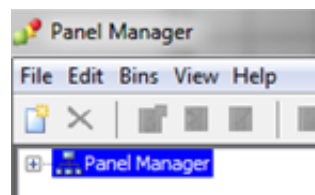
Before you use GeneMapper™ ID-X Software to analyze data for the first time, you must do the following:

**Set up GeneMapper™ ID-X Software**

- Check panel, bin, and stutter file versions on your computer**
- (If needed) Download newer versions of panel, bin, and stutter files**
- (If needed) Import panels, bins, and marker stutter**
- (Optional) Define custom table or plot settings**

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

1. Start the GeneMapper™ ID-X Software, then sign 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 the **Panels** folder, then check the version of panel, bin, and stutter files installed.



GeneMapper™ *ID-X* Software v1.7.x contains the latest panel, bin, and stutter files for the STR kits.

- If the latest files are not installed on your copy of the GeneMapper™ *ID-X* Software, proceed to “(If needed) Download newer versions of panel, bin, and stutter files” on page 32.
- If the latest files are already installed on your copy of the GeneMapper™ *ID-X* Software, skip to “Create an analysis method” on page 37.

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

1. Go to [www.thermofisher.com/GMIDXsoftware](http://www.thermofisher.com/GMIDXsoftware).  
The page provides a list of kit-specific analysis files. The analysis files for each kit can be downloaded in a single ZIP file.
2. If the analysis file versions listed for your kit are newer than the versions on your computer, download the ZIP file.

---

**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.

### **(If needed) Import panels, bins, and marker stutter**

Import the latest panel, bin set, and marker stutter from the website into the GeneMapper™ *ID-X* Software database.

---

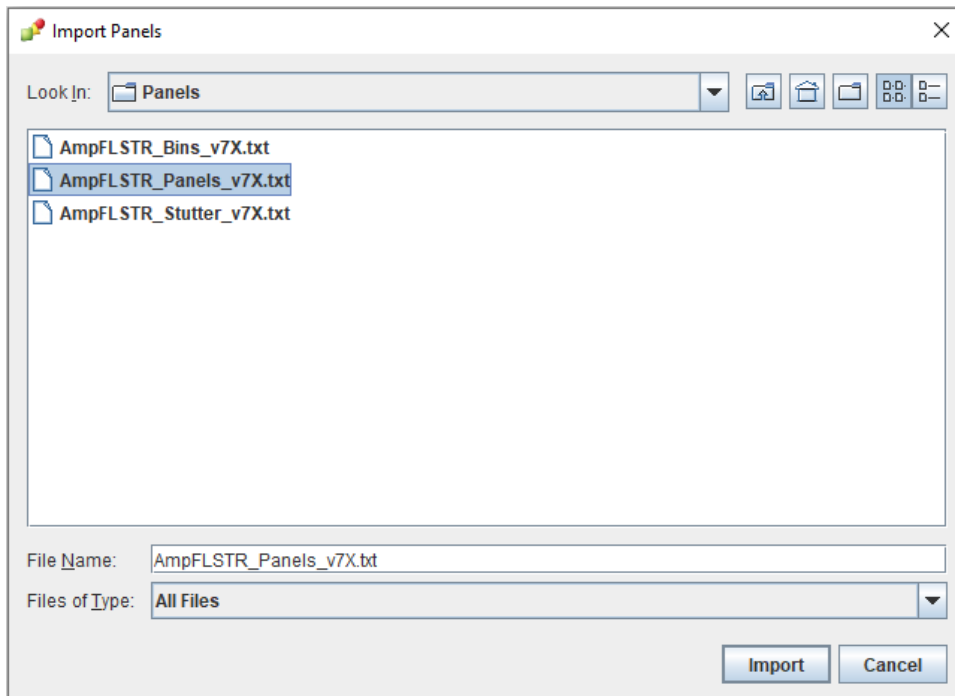
**Note:** The file names specified in this procedure are examples only. The files that you import may have different file names.

---

1. Start the GeneMapper™ *ID-X* Software, then sign in with the appropriate user name and password.
2. Select **Tools ▶ Panel Manager**.
3. Open the folder that contains 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 the analysis files folder that you unzipped in “(If needed) Download newer versions of panel, bin, and stutter files” on page 32.

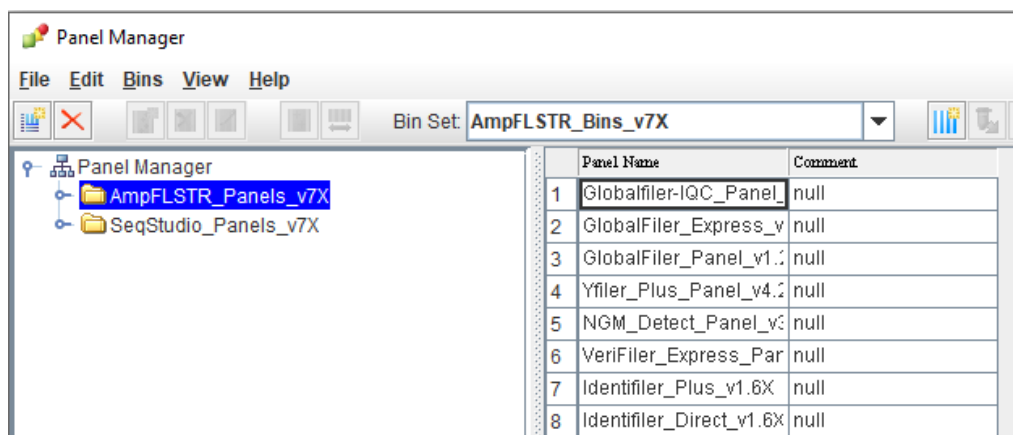


4. Select the panels TXT file for your kit, then click **Import**.



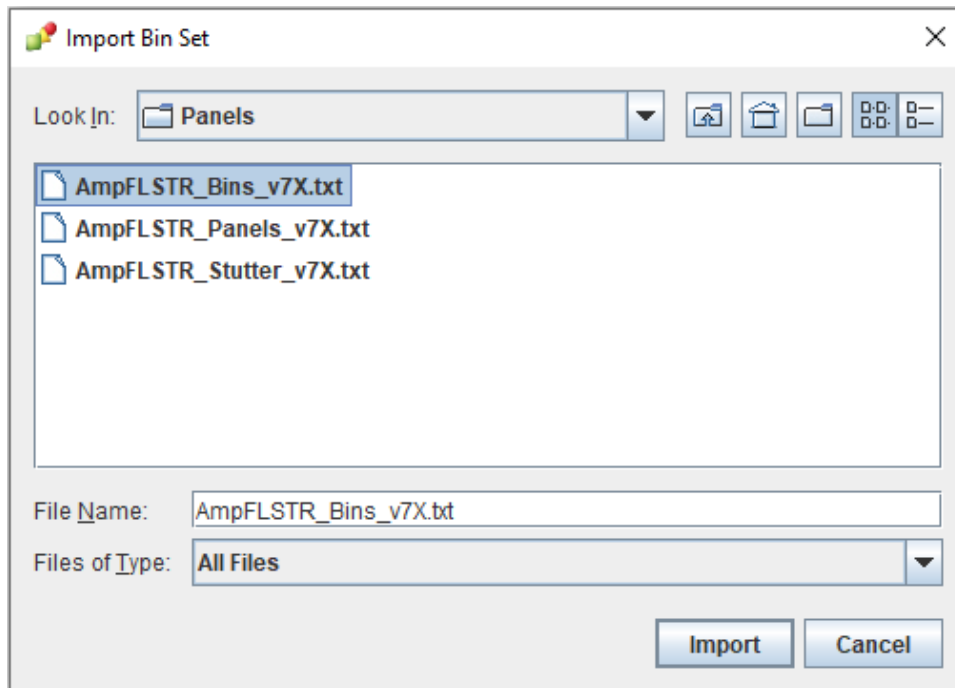
Importing the panels TXT file creates a new folder in the navigation pane of the **Panel Manager**. This folder contains the panels and associated markers.

5. Import the bins file.
  - a. In the navigation pane, select the panel folder created in step 4.



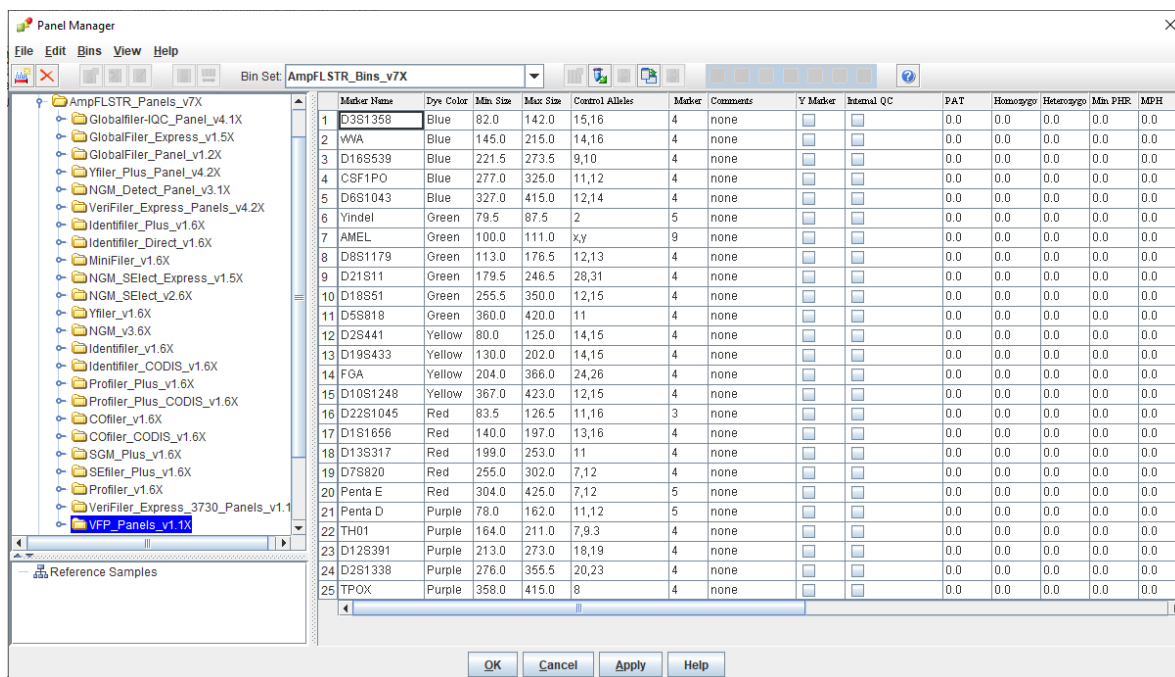
- b. Select **File** ► **Import Bin Set** to open the **Import Bin Set** dialog box.
- c. Navigate to the analysis files folder for your kit (from step 3).

- d. Select the bins TXT file for your kit, then click **Import**.

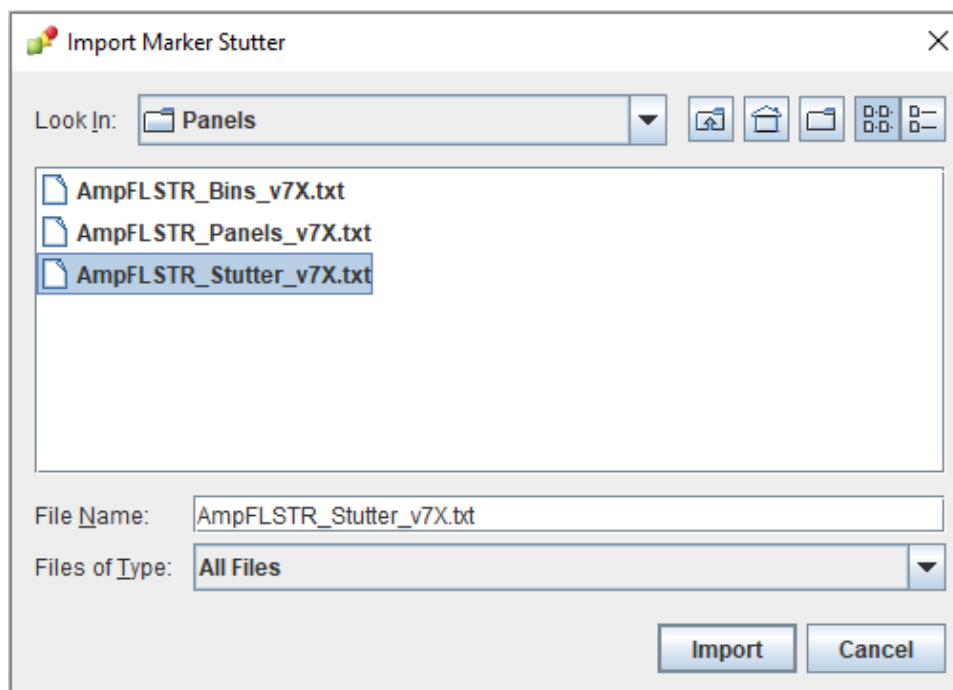


Importing the bins TXT file associates the bin set with the panels imported in step 4.

- 6. (Optional) View marker and panel information: In the navigation pane, select the panel folder for your kit. The panel information is displayed in the right pane and the markers are displayed below it.



7. Import the stutter file.
  - a. In the navigation pane, select the panel folder for your kit.
  - b. Select **File ▶ Import Marker Stutter** to open the **Import Marker Stutter** dialog box.
  - c. Navigate to the analysis files folder for your kit (from step 3).
  - d. Select the stutter TXT file for your kit, then click **Import**.



Importing the stutter TXT file associates the marker stutter ratio with the bin set in the panel folder for your kit (step 4) and overwrites any existing stutter ratios associated with the panels and bins in that folder.

8. (Optional) View the imported marker stutters.
  - a. In the navigation pane, click the panel folder for your kit to expand it.  
The markers are displayed in the navigation pane.

- b. Double-click a marker, then select the **Stutter Ratio & Distance** view for the marker in the right pane.

**Note:** The allele-specific stutter fields shown in the image are not implemented in GeneMapper™ ID-X Software v1.6 and earlier.

Panel Manager

File Edit Bins View Help

Bin Set: AmpFLSTR\_Bins\_v7X

Please enter the stutter filter(s) for D3S1358 marker here. If left blank, the global stutter filter will be applied.

**Marker Level Minus Stutter**

	Ratio	From Distance	To Distance
1	0.1261	3.25	4.75
2			
3			
4			

**Marker Level Plus Stutter**

	Ratio	From Distance	To Distance
1	0.0374	3.25	4.75
2			
3			
4			

**Allele-Specific Minus Stutter**

	Ratio	From Distance	To Distance	Allele
1				
2				
3				
4				

**Allele-Specific Plus Stutter**

	Ratio	From Distance	To Distance	Allele
1				
2				
3				
4				

New Edit Delete

New Edit Delete

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 GeneMapper™ ID-X Software database.

### (Optional) Define custom table or plot settings

Default views for table and plot settings are provided with the software.

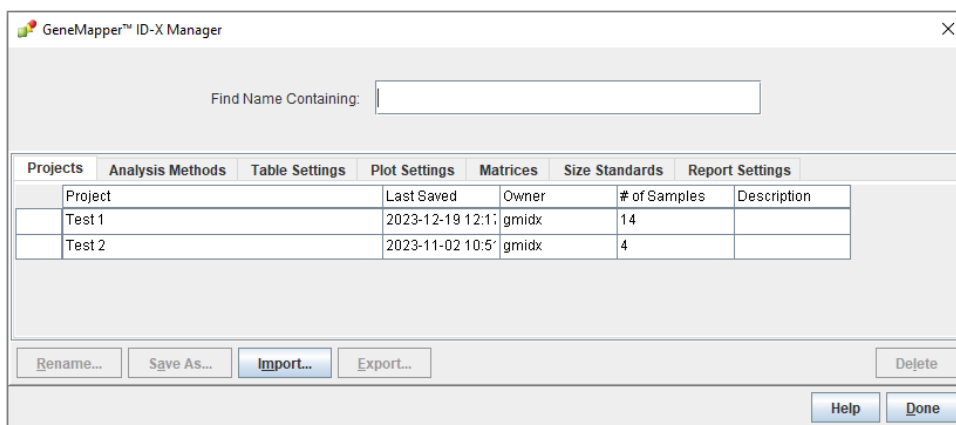
For information on defining custom views, see *GeneMapper™ ID-X Software v1.5 Getting Started Guide — Basic Features*.

## Create an analysis method

### Create an analysis method

**IMPORTANT!** Because analysis methods are version-specific, you need to create an analysis method for each version of the software. For example, an analysis method that is created in GeneMapper™ ID-X Software v1.6 is not compatible with analysis methods that are created in v1.5 or v1.7.x.

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 as described in the following pages.

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

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

## Enter Analysis Method settings

### Enter General tab settings

1. Enter an analysis method name.
2. Select the security group appropriate for your software configuration.
3. *(Optional)* Enter a description and an instrument.

The screenshot shows the 'Analysis Method Editor' dialog box with the 'General' tab selected. The dialog has a title bar with a close button (X) and a tabbed interface with four tabs: 'General', 'Allele', 'Peak Detector', and 'SQ & GQ Settings'. The 'General' tab is active and contains the following fields:

- Analysis Method Description** (Section Header)
- Name:** My\_Analysis\_Method
- Security Group:** GeneMapper ID-X Security Group (dropdown menu)
- Description:** (empty text area with scrollbars)
- Instrument:** (empty text field)
- Analysis Type:** HID

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

Figure 6 General tab settings

## Enter Allele tab settings

**IMPORTANT!** Perform internal validation studies to determine the appropriate settings for your laboratory.

1. Select the appropriate bin set.
2. (Optional) Select stutter options.

Option	Action	Additional information
<b>Use marker-specific stutter ratio and distance if available</b>	Select or deselect the checkbox, as needed.	To apply the stutter ratios that are contained in the Panel Manager, select the checkbox.
<b>Use allele-specific stutter ratios and distances if available</b> The checkbox is available only for GeneMapper™ ID-X Software v1.7 or later.	Select or deselect the checkbox, as needed.	To use allele-specific stutter filtering, select the checkbox.
<b>Consider additive stutters (forward and back)</b> The checkbox is available only for GeneMapper™ ID-X Software v1.7 or later.	Select or deselect the checkbox, as needed.	To take additive stutter into consideration, select the checkbox.

**Note:** For more information on the GeneMapper™ ID-X Software v1.7 options, see the *GeneMapper™ ID-X Software v1.7 New Features and Software Verification and Validation User Bulletin* (Pub. No. [MAN0029209](#)).

3. In the **Marker Repeat Type** pane, enter values for the Tri, Tetra, Penta, and Hexa loci.

**Note:** For paternity and database applications: In the **Global Cut-off Value** field, we recommended using a cut-off value of 20% for the Tri, Tetra, and Penta loci.

4. Enter the appropriate filter settings.

**Analysis Method Editor** [X]

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

Bin Set: **AmpFLSTR\_Bins\_v7X** ▼

Use marker-specific stutter ratio and distance if available  
 Use allele-specific stutter ratios and distances if available.  
 Consider additive stutters (forward and back).

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

Amelogenin Cutoff

Figure 7 Allele tab settings



## Enter Peak Detector tab settings

Enter or select the appropriate values.

Option	Action	Additional information
<b>Use marker-specific thresholds (if available)</b> The checkbox is available only for GeneMapper™ ID-X Software v1.7 or later.	Select or deselect the checkbox, as needed.	To use the marker-specific thresholds defined in the Panel Manager, select the checkbox.
<b>Ranges</b>	<b>Analysis</b> —Select <b>Full Range</b> from the dropdown list.	<i>(if needed)</i> The analysis range can be narrowed.
	<b>Sizing</b> —Select <b>All Sizes</b> from the dropdown list.	—
<b>Smoothing and Baselineing</b>	<b>Smoothing</b> —Select <b>Light</b> .	The Verifiler™ Plus kit was validated with the values listed. If your laboratory uses an unvalidated polymer, you may need to adjust these values.
	<b>Baseline Window</b> —Enter <b>33</b> pts.	
<b>Size Calling Method</b>	Select <b>Local Southern Method</b> or <b>3rd Order Least Squares</b> .	The Verifiler™ Plus kit was validated using the Local Southern Method and 3rd Order Least Squares. Do not select another method unless you perform internal validation studies to determine the appropriate method for your laboratory.
<b>Peak Detection</b>	<b>Peak Amplitude Thresholds</b> —User-defined.	The default value is 50 RFU for all dyes. Perform internal validation studies to determine the appropriate peak amplitude thresholds for your laboratory.
	<b>Min. Peak Half Width</b> —Enter <b>2</b> pts.	The Verifiler™ Plus kit was validated with the values listed. Do not enter other values unless you perform internal validation studies to determine the appropriate values for your laboratory.
	<b>Polynomial Degree</b> —Enter <b>3</b> .	
	<b>Peak Window Size</b> —Enter <b>11</b> pts.	
<b>Slope Threshold</b>	<b>Peak Start</b> —Enter <b>0.0</b> .	
	<b>Peak End</b> —Enter <b>0.0</b> .	

(continued)

Option	Action	Additional information
<p><b>Use Normalization, if applicable</b></p> <p>The checkbox is available for use with data run on the following instruments:</p> <ul style="list-style-type: none"> <li>• SeqStudio™ Flex Series Genetic Analyzer for Human Identification</li> <li>• 3500 Series Genetic Analyzer for Human Identification</li> </ul>	<p>Select or deselect the checkbox, as needed.</p>	<p>To apply size standard normalization data to the analysis, select the checkbox.</p> <p>The size standard normalization data are collected on the capillary electrophoresis instrument. To see if normalization data have been collected for a specific data file, see <b>SS Normalization Factor</b> in the GeneMapper™ ID-X Software.</p>

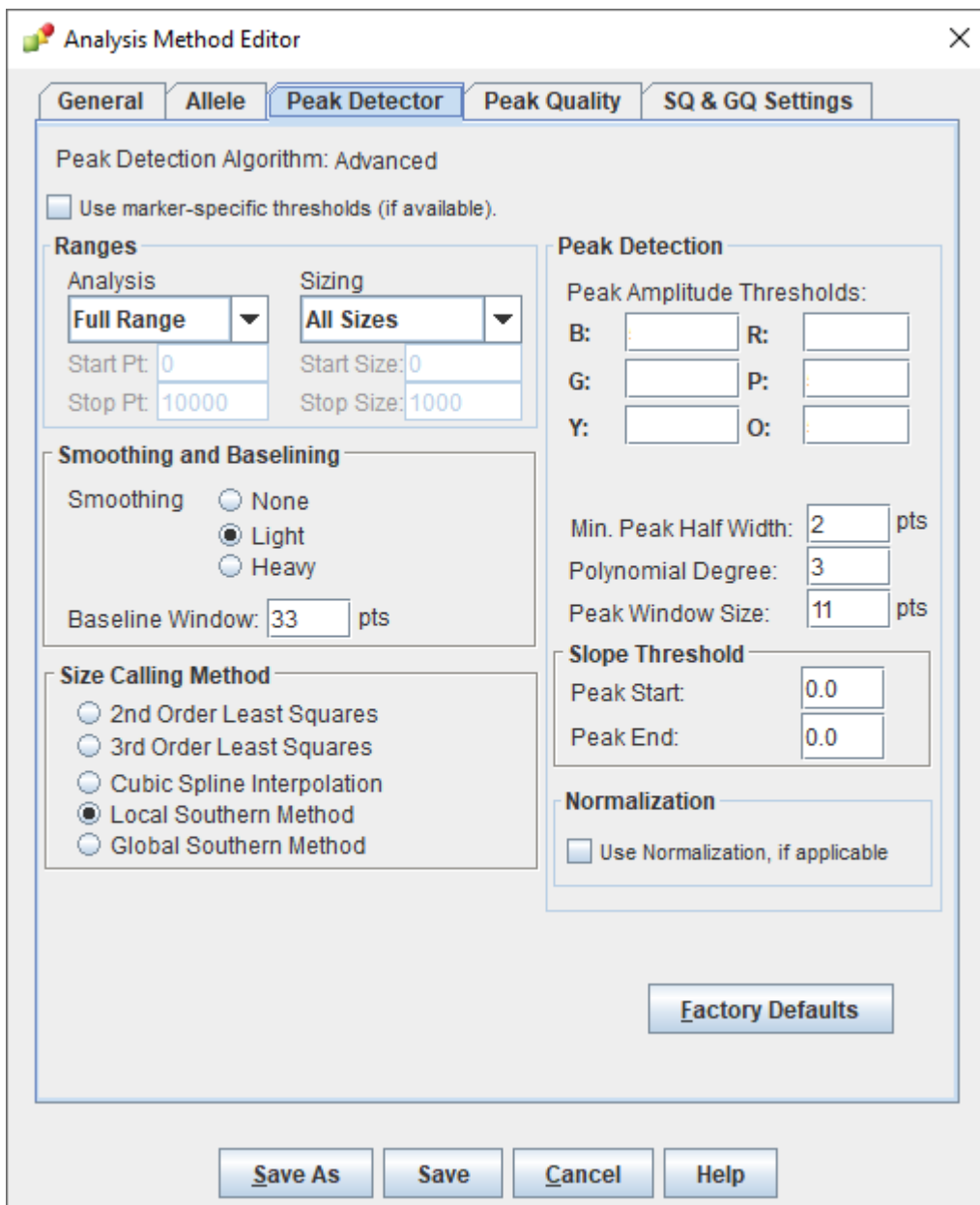


Figure 8 Peak Detector tab settings

### Enter Peak Quality tab settings

1. Perform 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.
2. For the remaining fields, enter the values shown in Figure 9.

**Note:** The **Pull-Up Ratio (PU)** settings shown in the figure are implemented only in GeneMapper™ ID-X Software v1.7 or later. For more information on these settings, see the *GeneMapper™*

ID-X Software v1.7 New Features and Software Verification and Validation User Bulletin  
(Pub. No. [MAN0029209](#)).

The screenshot shows the 'Analysis Method Editor' window with the 'Peak Quality' tab selected. The settings are as follows:

Section	Parameter	Value
Min/Max Peak Height (LPH/MPH)	Homozygous min peak height	[Empty]
	Heterozygous min peak height	[Empty]
	Max Peak Height (MPH)	[Empty]
Peak Height Ratio (PHR)	Min peak height ratio	[Empty]
Broad Peak (BD)	Max peak width (basepairs)	1.5
Allele Number (AN)	Max expected alleles:	
	For autosomal markers & AMEL	2
	For Y markers	1
Allelic Ladder Spike	Spike Detection	Enable
	Cut-off value	0.2
Sample Spike Detection	Spike Detection	Enable
Pull-Up Ratio (PU)	Enable pull-up detection.	<input type="checkbox"/>
	Label pull-up	<input checked="" type="radio"/>
	Remove pull-up peaks	<input type="radio"/>
	Max pull-up ratio	0.05
Pull-up offset (data points)		2

Buttons at the bottom: Save As, Save, Cancel, Help.

Figure 9 Peak Quality tab settings

### Enter SQ & GQ tab settings

Enter the appropriate values.

**IMPORTANT!** The software default values are shown in Figure 10. We used the software default values during developmental validation. We recommend that you perform internal validation studies to determine the appropriate values for your laboratory.

**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.

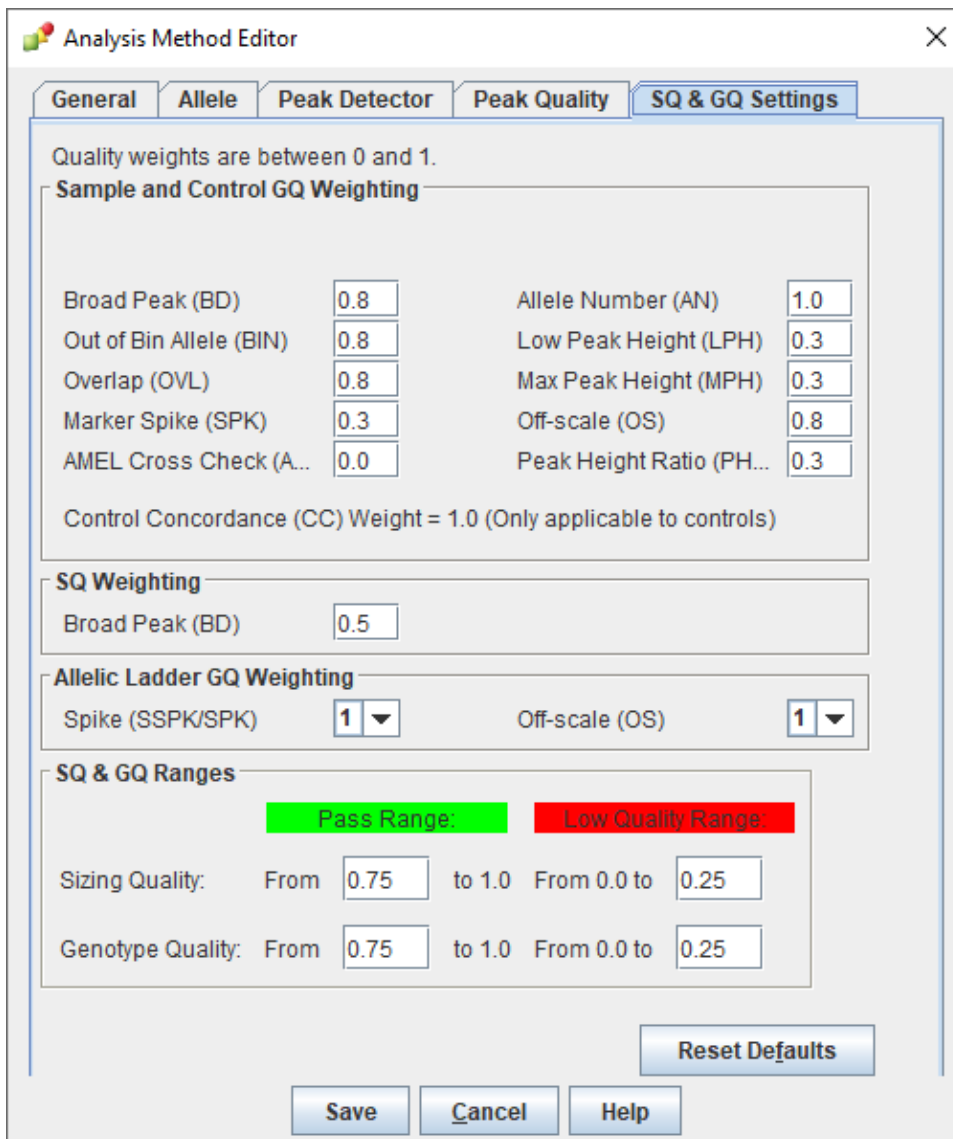


Figure 10 SQ & GQ tab settings

## (If needed) Create a size standard definition file

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

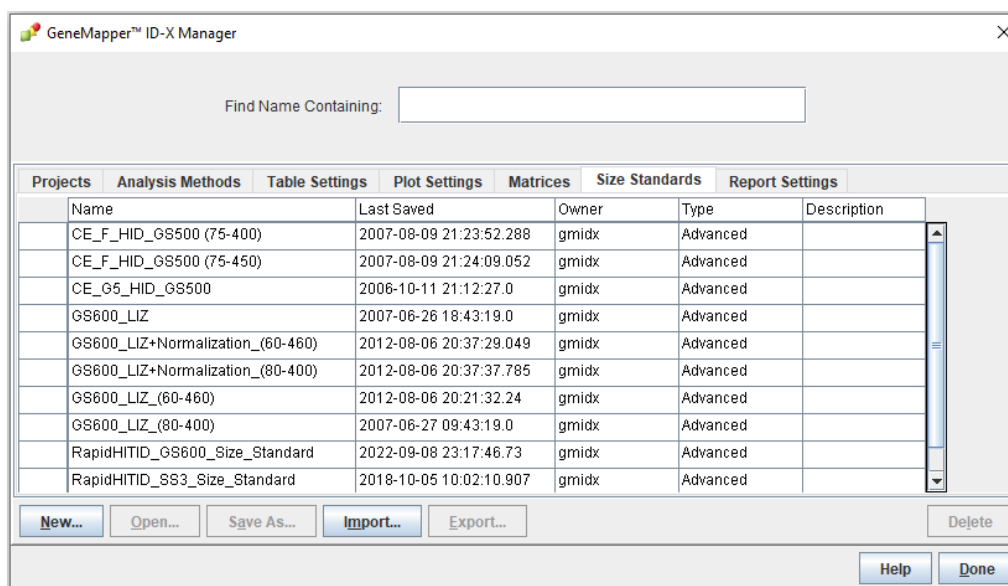
## About the GS600\_LIZ\_(60–460) size standard definition file

The GS600\_LIZ\_(60–460) size standard definition file that is provided with GeneMapper™ ID-X Software and used with the Local Southern Method (size calling method) contains the following peaks: 60, 80, 100, 114, 120, 140, 160, 180, 200, 214, 220, 240, 250, 260, 280, 300, 314, 320, 340, 360, 380, 400, 414, 420, 440, and 460.

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 46.

## 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.
  - 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. (If needed) Click **Insert** to add rows or click **Delete** to remove rows.

**Size Standard Editor**

**Edit**

**Size Standard Description**

Name: My\_Size\_Standard

Security Group: GeneMapper ID-X Security Group

Description:

Size Standard Dye: Orange

**Size Standard Table**

	Size in Basepairs
1	0.0
2	0.0
3	0.0
4	0.0
5	0.0
6	0.0
7	0.0

Insert Delete

OK Cancel Help

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

1. 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.

Option	Action
<b>Sample Type</b>	Select the sample type for each sample, control, and allelic ladder in the project.
<b>Analysis Method</b>	Select the analysis method that you created in “Create an analysis method” on page 37.
<b>Panel</b>	Select the current kit panel. If needed, see “Check panel, bin, and stutter file versions on your computer” on page 31.
<b>Size Standard</b>	Select the <b>GS600_LIZ_(60-460) size standard definition</b> , or select another validated size standard definition, as described in “(If needed) Create a size standard definition file” on page 45.

3. Click **Analyze**.
4. In the **Save Project** dialog box, enter a name for the project, then click **OK** to start analysis.
  - The status bar displays the progress of analysis.
  - 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™ ID-X Software

See “Related documentation” on page 133 for a list of available documents.





# Assess the PCR reaction with the Internal Quality Control System

■ Overview of the Internal Quality Control system .....	49
■ Evaluate the PCR reaction .....	49

## Overview of the Internal Quality Control system

The Internal Quality Control (IQC) System is a tool that can help you evaluate the PCR reaction and, along with the STR marker data, infer possible sample degradation or inhibition.

The primers for the two IQC markers, IQCS and IQCL, amplify synthetic DNA targets that are included in the primer set. IQCS is a low molecular weight amplicon, with mobility of 70 bp. IQCL is a higher molecular weight amplicon, with mobility of 451 bp.

---

**Note:** The IQC markers enable *qualitative* insight into the sample amplification. By default, the IQC markers are not genotyped during analysis in the GeneMapper™ *ID-X* Software. However, if needed, the Verifiler™ Plus Allelic Ladder contains two peaks each for the IQCS and IQCL markers (designated 1 and 2) to detect the IQC sequences. Sample reactions contain one peak each for IQCS and IQCL, which corresponds to the larger peak (peak 2) in each IQC marker.

---

The IQC system allows 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 might be present in the PCR reaction, or if the PCR reaction conditions are not optimal, by evaluating the peak heights of IQCS and IQCL.

## Evaluate the PCR reaction

To evaluate the PCR performance of the samples, review the peak heights of the IQCS and IQCL. Under ideal PCR conditions, the peak heights of the IQCS and IQCL should be >2,000 RFU using standard injection protocols on the 3500/3500xL Genetic Analyzer (see Table 5). Under suboptimal PCR conditions (for example, moderate inhibition), the height of the IQCL is substantially reduced. Under extremely high inhibition, even the small IQC peak heights are substantially reduced. Note that when high inputs of DNA are amplified (>2 ng) some suppression of the IQCL peak may also be seen.

For outcome scenarios, see Table 5.

**Note:** Electropherograms from runs on a 3500xL Genetic Analyzer using standard injection protocols are shown in Figure 11 through Figure 16.

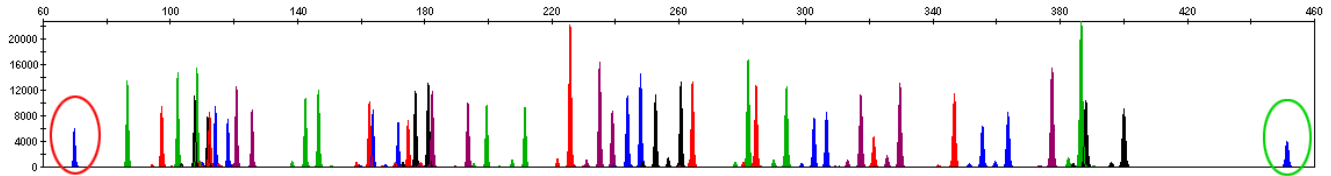
**Table 5 IQC peak interpretation**

Sample DNA profile	IQC peaks	IQC interpretation	Recommended action for re-amplification <sup>[1]</sup>
		Typically indicates	
Balanced	Both IQC peaks are >2,000 RFU	PCR performance within the optimal range—no sample degradation or inhibition (see Figure 11)	Re-amplification should not be required. On a case-by-case basis, you can evaluate the profile and determine if re-amplification is required.
Balanced	The IQCS peak height is >2,000 RFU, while the IQCL peak height is significantly decreased	Mild inhibition (see Figure 12)	Evaluate the profile and determine if additional input would recover alleles without saturating the system.
Ski-slope	The IQCS peak height is near 2,000 RFU or greater, while the IQCL peak is not present	High inhibition (see Figure 13)	Perform a dilution or dilutions of the sample to minimize the inhibitor while still allowing for enough DNA to achieve amplification (for example, 10–20% dilution).
Ski-slope	The IQCS peak height is significantly decreased, while the IQCL peak is not present	Very high inhibition (see Figure 14)	Perform a dilution or dilutions of the sample to minimize the inhibitor while still allowing for enough DNA to achieve amplification (for example, 20–60% dilution).
Ski-slope	Both IQC peaks are >2,000 RFU	Degraded sample DNA (see Figure 15)	Evaluate the profile and determine if additional input would recover alleles without saturating the system.
No peaks	Both IQC peaks are >2,000 RFU	No DNA or very little sample DNA (see Figure 16)	Check the quantification and normalization calculations. If you have sufficient DNA to generate a profile, increase the amount of DNA in the amplification.
No peaks	No peaks	PCR failure or ultra-high inhibition	Re-amplify the sample at the appropriate target to determine if PCR failure occurred. If DNA is still not detected, perform a dilution or dilutions of the sample to minimize the inhibitor while still allowing for enough DNA to achieve amplification (for example, 20–60% dilution).

<sup>[1]</sup> The decision to re-amplify the sample should be based on individual laboratory protocols that identify how much information is sufficient for reporting.

## Balanced profile

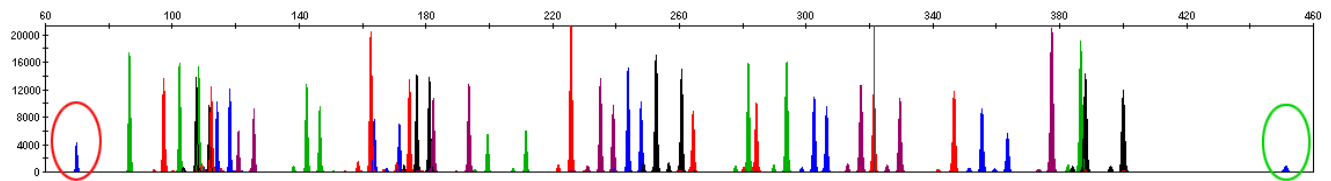
In this example of a balanced profile, the IQC peaks and the DNA profile peaks are balanced, which indicates that PCR has occurred optimally.



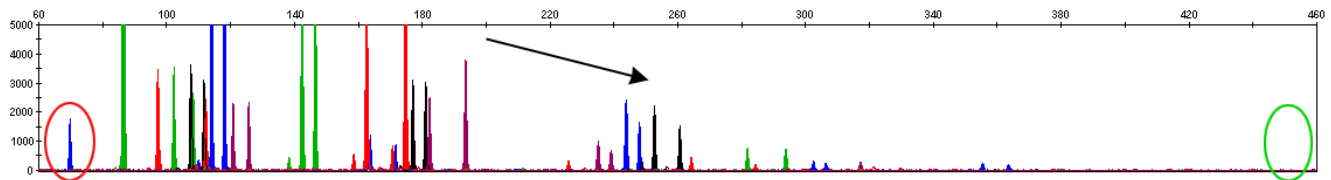
**Figure 11** Combined dyes electropherogram for the Verifiler™ Plus kit with 0.5 ng of DNA (Y-axis scale 0–22,000 RFU) (updated formulation). The red circle highlights the FAM™ dye channel IQCS peak and the green circle highlights the FAM™ dye channel IQCL peak.

## Ski slope profile with decreased IQCL peak height

A significantly lower IQCL peak height is shown in Figure 12. This indicates that the PCR reaction has been compromised by inhibition. The complete absence of an IQCL peak, indicating a high level of inhibition, is shown in Figure 13.



**Figure 12** Combined dyes electropherogram for the Verifiler™ Plus kit in the presence of 150 ng/μL of humic acid (Y-axis scale 0–22,000 RFU) (updated formulation). The red circle highlights the IQCS peak (a PCR inhibitor) and the green circle highlights the IQCL peak. While overall peak heights are reduced relative to the uninhibited control sample, only a subtle ski-slope effect is observed.



**Figure 13** Combined dyes electropherogram for the Verifiler™ Plus kit in the presence of 350 ng/μL of humic acid (Y-axis scale 0–5,000 RFU) (updated formulation). The red circle highlights the IQCS peak; the IQCL peak is absent (green box). The arrow indicates the ski slope peak pattern observed in the DNA profile.

## Ski slope profile with IQC peaks

The presence of both IQCS and IQCL peaks >2,000 RFU indicates that PCR has occurred optimally.

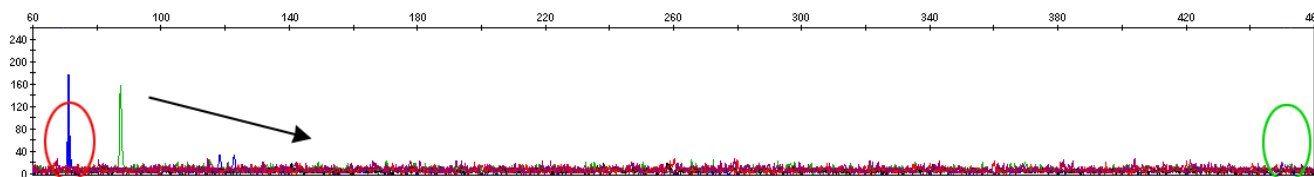


Figure 14 Combined dyes electropherogram for the Verifiler™ Plus kit in the presence of 450 ng/μL of humic acid (updated formulation). The red circle highlights the IQCS peak; the IQCL peak is absent (green circle). The arrow indicates the ski slope peak pattern observed in the DNA profile. The Y-axis is scaled to 250 RFU to make extremely low peaks visible.

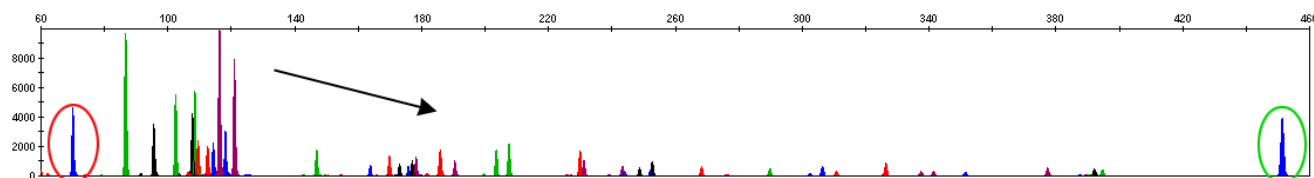


Figure 15 Combined dyes electropherogram from degraded DNA (Y-axis scale 0–10,000 RFU) (updated formulation). The red circle highlights the IQCS peak and the green circle highlights the IQCL peak. The arrow indicates the ski slope pattern observed in the DNA profile.

## No sample peaks with IQC peaks

There are no DNA profile peaks in Figure 16. However, the presence of both the IQCS and IQCL peaks indicates that normal amplification has occurred in the PCR.

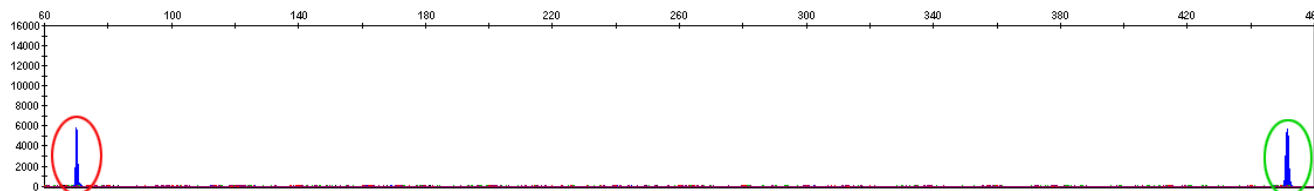


Figure 16 Combined dyes electropherogram that shows IQCS and IQCL peaks with 0 ng of DNA (Y-axis scale 0–16,000 RFU) (updated formulation). The red circle highlights the IQCS peak and the green circle highlights the IQCL peak.



# Experiments and results

- Importance of validation ..... 53
- Experiment conditions ..... 53
- Laboratory requirements for internal validation ..... 54
- Developmental validation ..... 54
- Accuracy, precision, and reproducibility ..... 57
- Extra peaks in the electropherogram ..... 76
- Characterization of loci ..... 85
- Species specificity ..... 87
- Sensitivity ..... 89
- Stability ..... 90
- Mixture studies ..... 94
- Population data ..... 98

## 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 (SWGDM, 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.

## Repeated validation experiments using the updated primer set formulation

In 2020, minor modifications were made to the Verifiler™ Plus PCR Amplification Kit primer set to decrease the incidence of non-specific PCR artifacts that were observed in certain forensic samples with high microbial DNA load. Accordingly, we used the updated primer set formulation to repeat the validation experiments that might be impacted by the primer set changes, as follows:

- Developmental validation—“PCR components” on page 55
- Extra peaks—“Dye artifact observation” on page 83
- Species Specificity—“Nonhuman study observation” on page 87
- Sensitivity—“Sensitivity observation” on page 89
- Stability—“Degraded DNA” on page 91
- Stability—“Effect of inhibitors” on page 91
- Mixture Studies—“Limit of detection of the minor component” on page 96

Because most kit components (for example, the master mix) remain unchanged, we did not repeat the remaining validation experiments.

---

**Note:** In this chapter, "updated formulation" refers to the 2020 updated primer set formulation.

---

## 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

Developmental validation studies were performed using the ProFlex™ 96-well PCR System, Veriti™ 96-Well Thermal Cycler, and the GeneAmp™ PCR System 9700 96-Well thermal cycler according to the protocol described in the Perform PCR chapter.

### SWGAM 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.” (SWGAM, December 2016)

### SWGAM 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.” (SWGAM, December 2016)

## SWGAM guideline 3.9.6

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

## PCR components

We examined the concentration of each component in the kit. We established that the concentration of each component was within the range where data indicated that the amplification met the required performance criteria for specificity, sensitivity, and reproducibility.

For example, 0.5 ng of DNA Control 007 was amplified in the presence of varying concentrations of magnesium sulfate, and the results were analyzed on a 3500xL Genetic Analyzer (Figure 17). The performance of the multiplex is most robust within  $\pm 20\%$  of the optimal magnesium sulfate concentration.

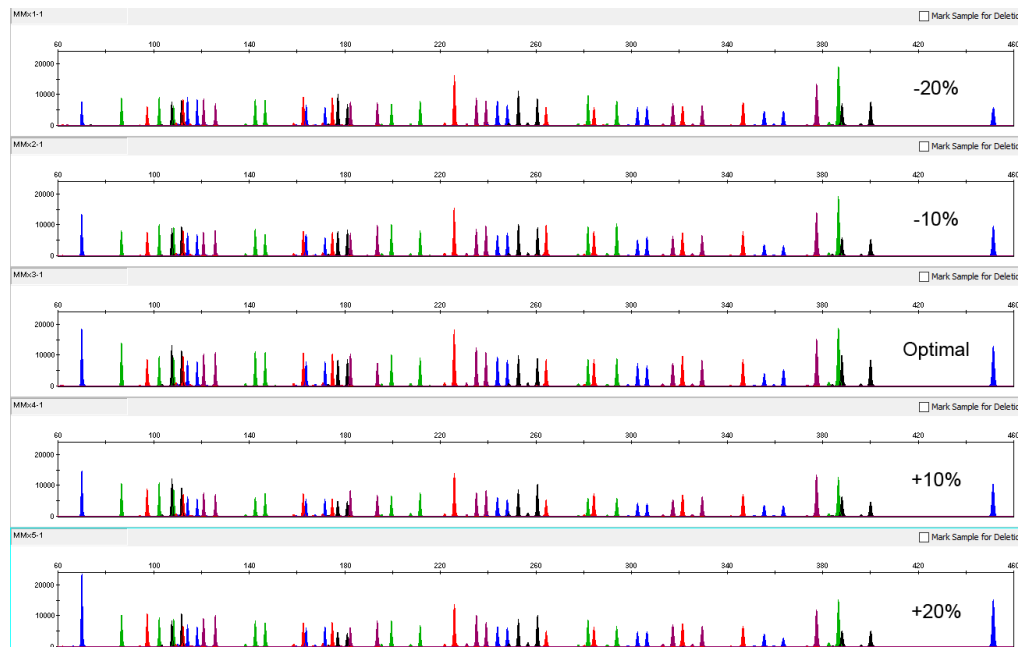


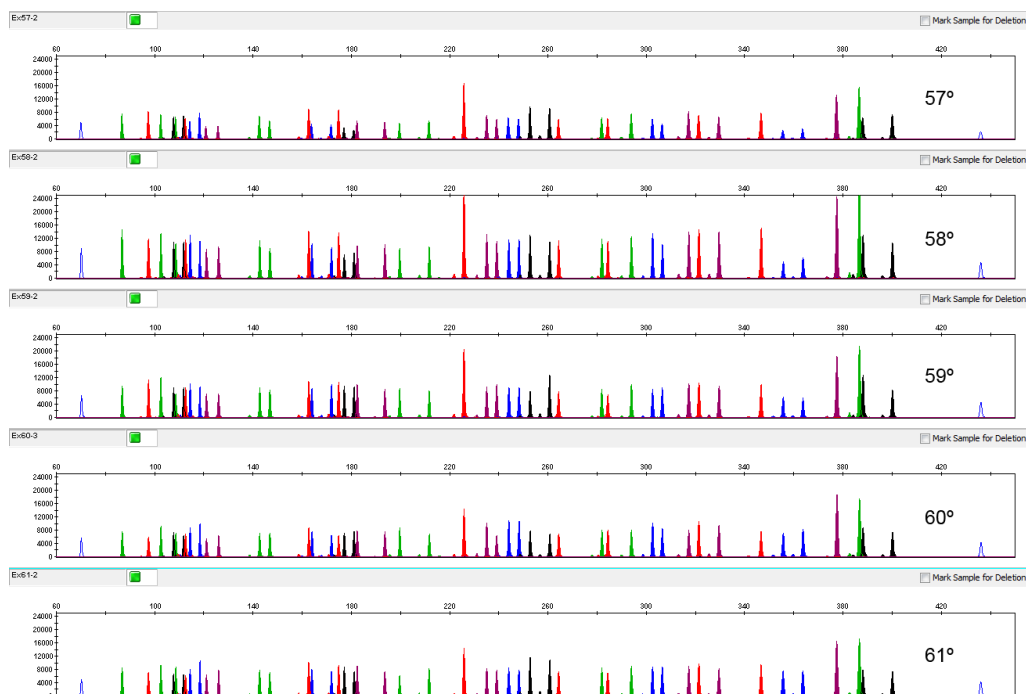
Figure 17 DNA Control 007 (0.5 ng) amplified with the Verifiler™ Plus kit in the presence of varying concentrations of magnesium sulfate and analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0–25,000 RFU) (updated formulation)

## Thermal cycling temperatures

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

For example, annealing/extension temperatures of 57, 58, 59, 60, and 61°C were tested using a ProFlex™ PCR System (Figure 18). The PCR products were analyzed using a 3500xL Genetic Analyzer.

Robust profiles were obtained between 57–61°C. 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.



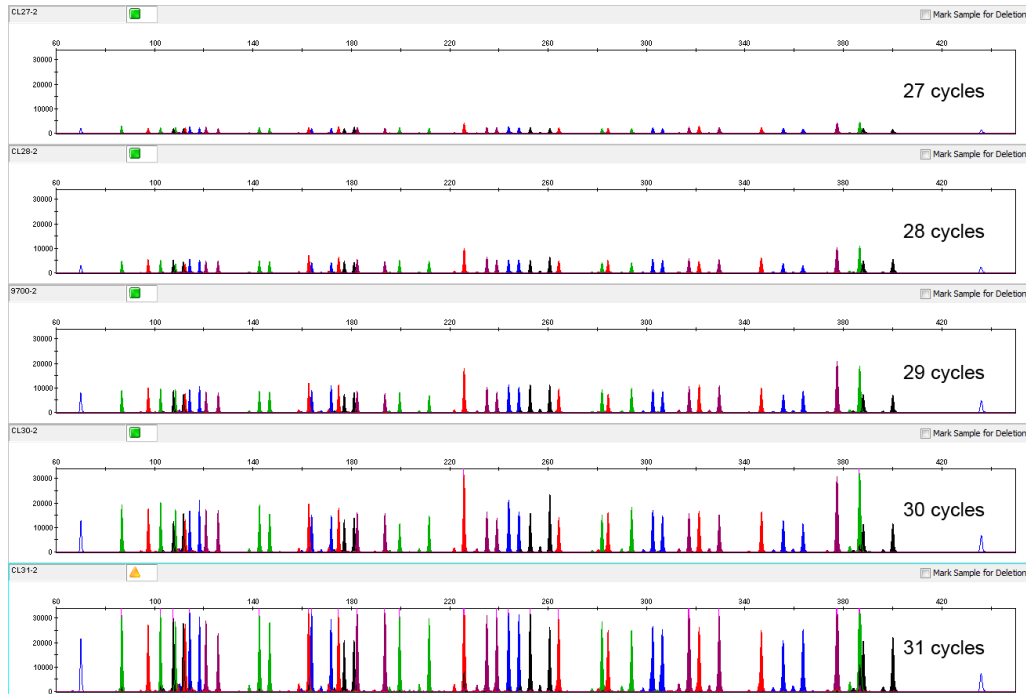
**Figure 18** Verifiler™ Plus kit electropherograms obtained from amplification of 0.5 ng of DNA Control 007 at annealing temperatures of 57, 58, 59, 60, and 61°C, analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0–25,000 RFU)

## PCR cycle number

Reactions were amplified for 27, 28, 29, 30, and 31 cycles on the ProFlex™ PCR System 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 (27–31) and off-scale data were collected for several allele peaks at 31 cycles (Figure 19).

Additional experiments were performed not only to optimize for peak heights, peak balance, sensitivity, and so on, but also to minimize the formation of non-target PCR artifacts. The optimized end result is a "touchdown" PCR protocol in which the first two cycles are performed at a slightly higher, more stringent annealing temperature (62°C), followed by 27 cycles with an annealing temperature of 59°C.





**Figure 19** Representative Verifiler™ Plus kit profiles obtained from amplification of 0.5 ng of DNA Control 007 using 27, 28, 29, 30, and 31 cycles, analyzed on a 3500xL Genetic Analyzer (Y-axis scale 0–35,000 RFU).

## Accuracy, precision, and reproducibility

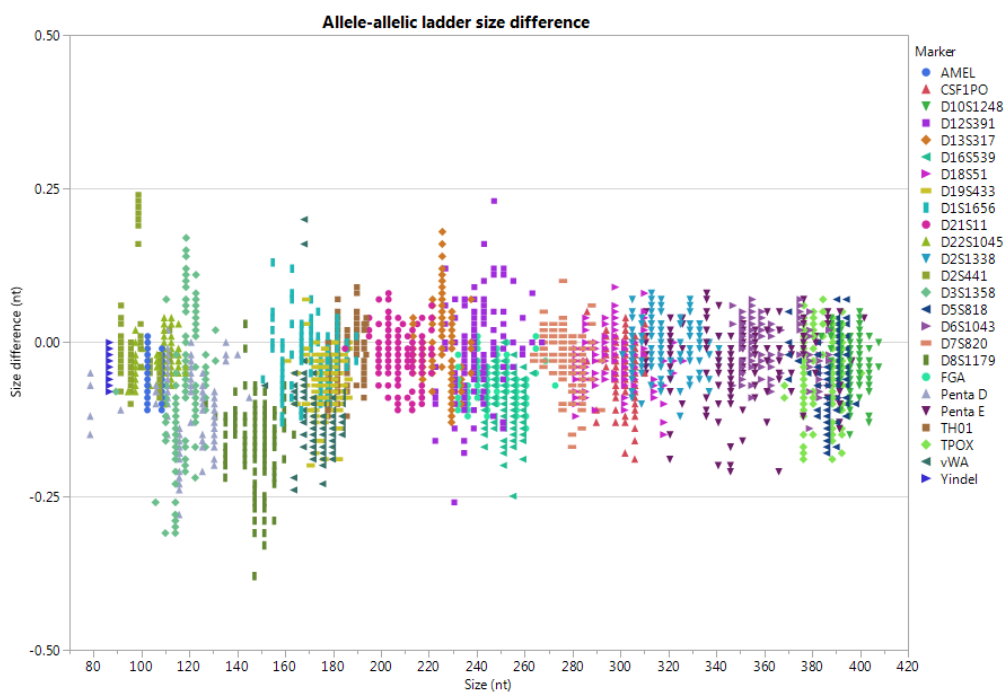
### SWGDM guideline 3.5

“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.” (SWGDM, December 2016)

### Accuracy observation

The size differences that were observed between sample alleles and the Verifiler™ Plus Allelic Ladder alleles on the 3130xL, 3500, and 3500xL instruments with POP-4™ Polymer are shown in Figure 20, Figure 21, and Figure 22. The X-axis represents the nominal base pair sizes for the allelic ladder. The horizontal lines parallel to the X-axis represent the  $\pm 0.25$ -bp windows. The Y-axis represents the deviation of each sample allele size from the corresponding allelic ladder allele size. All sample alleles are within  $\pm 0.5$  bp from a corresponding allele in the allelic ladder, regardless of the capillary electrophoresis platform.

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



**Figure 20** Allele size vs. allelic ladder sizing for 84 samples analyzed on a 3130xl Genetic Analyzer. Size and ladder sizing for the Verifiler™ Plus kit were calculated using the GeneScan™ 600 LIZ™ Size Standard v2.0.

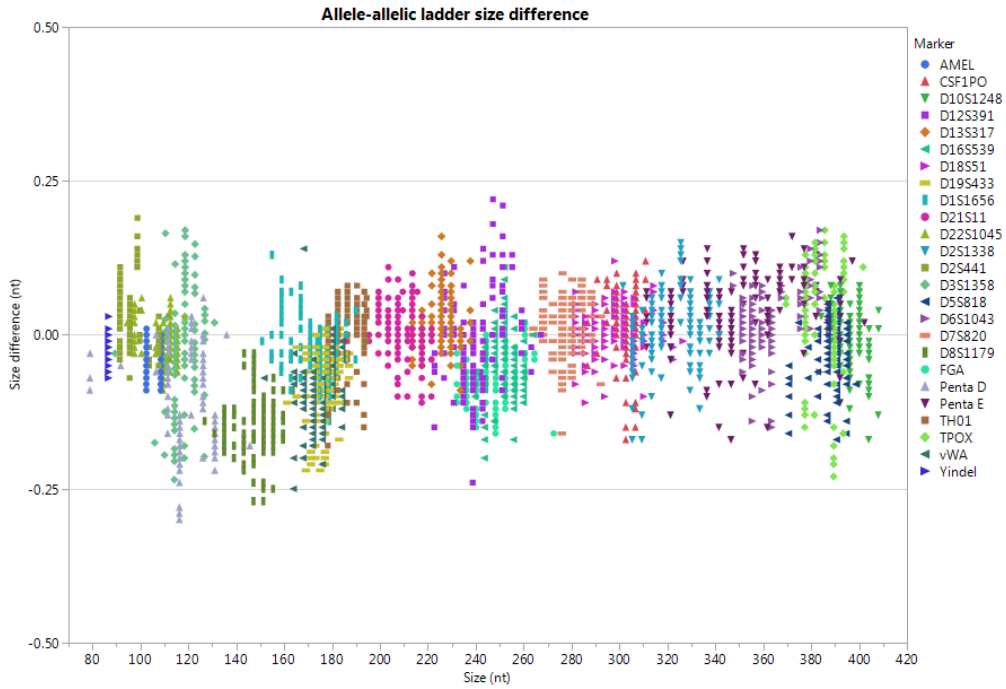


Figure 21 Allele size vs. allelic ladder sizing for 84 samples analyzed on a 3500 Genetic Analyzer. Size and ladder sizing for the Verifiler™ Plus kit were calculated using the GeneScan™ 600 LIZ™ Size Standard v2.0.

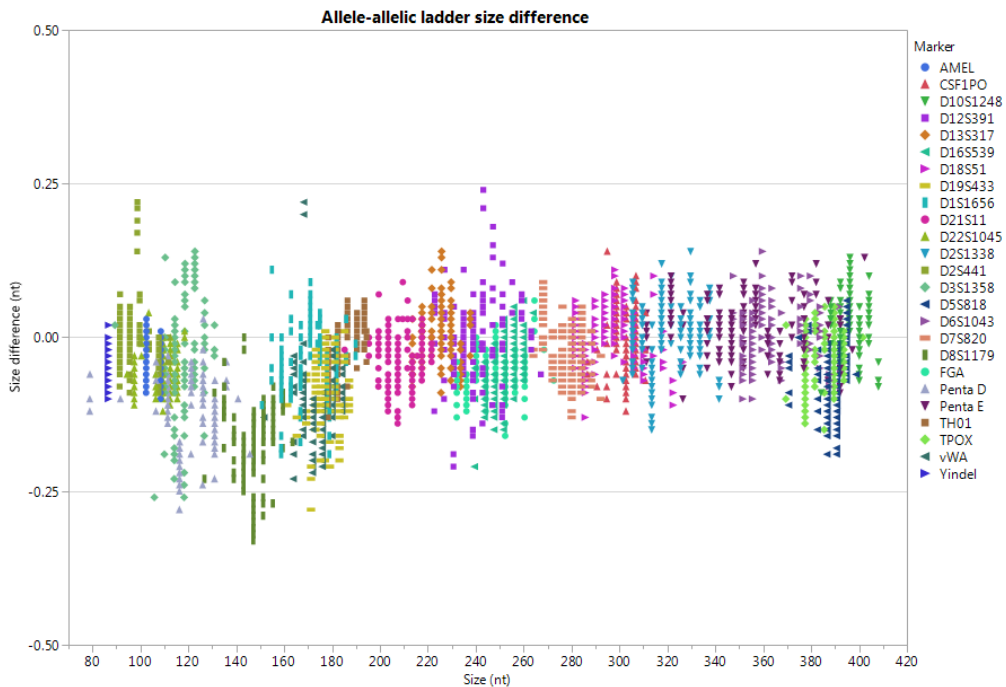


Figure 22 Allele size vs. allelic ladder sizing for 84 samples analyzed on a 3500xL Genetic Analyzer. Size and ladder sizing for the Verifiler™ Plus kit were calculated using the GeneScan™ 600 LIZ™ Size Standard v2.0.

## Precision and size window description

Sizing precision enables the determination of accurate and reliable genotypes. The recommended method for genotyping is to use a  $\pm 0.5$ -bp “window” around the size obtained for each allele in the allelic ladder. A  $\pm 0.5$ -bp window allows for the detection and correct assignment of alleles. Any sample allele that sizes outside the specified window could be one of the following:

- 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 electrophoresis instrument.

## Precision and size window observation

Typical precision results obtained from multiple runs of the Verifiler™ Plus Allelic Ladder using the GeneScan™ 600 LIZ™ Size Standard v2.0 are shown in Table 7. The results were obtained within a set of injections on a single capillary array. The number of repeated injections for each capillary electrophoresis platform is shown in Table 6.

**Table 6** Repeated injections for each CE platform

Genetic analyzer	Capillaries	Repeated injections	Sizing method
3130xl	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 7). 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  $\pm 0.5$ -bp window for a respective allelic ladder allele because of measurement error. The frequency of such an occurrence is lowest in detection systems with the smallest standard deviations in sizing. The tight clustering of allele sizes obtained, where the standard deviation in sizing is typically  $< 0.15$  bp, is illustrated in Figure 20 through Figure 22. The instance of a sample allele sizing outside the  $\pm 0.5$ -bp window because of measurement error is relatively rare when the standard deviation in sizing is approximately  $\leq 0.15$  bp (Smith, 1995).

For sample alleles that do not size within a  $\pm 0.5$ -bp window, the PCR product must be rerun to distinguish between a true off-ladder allele versus measurement error of a sample allele that corresponds to an allele in the allelic ladder. Repeat analysis, when necessary, provides an added level of confidence in the final allele assignment.

The GeneMapper™ *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 help ensure accurate genotyping, follow the guidelines in “Allelic ladder requirements for data analysis” on page 30 and use allelic ladders obtained from the same run as samples to analyze the samples.

**Note:** 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 7 Precision results of multiple runs of the Verifiler™ Plus Allelic Ladder

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
<b>D3S1358</b>						
9	89.28–89.33	0.013–0.046	89.48–89.52	0.006–0.039	89.37–89.44	0.010–0.045
10	93.45–93.52	0.012–0.039	93.67–93.70	0.005–0.046	93.55–93.63	0.013–0.039
11	97.62–97.69	0.012–0.042	97.83–97.89	0.006–0.042	97.75–97.79	0.008–0.054
12	101.78–101.86	0.018–0.040	102.00–102.06	0.010–0.040	101.91–101.96	0.010–0.043
13	105.95–105.99	0.014–0.041	106.13–106.20	0.010–0.040	106.04–106.11	0.008–0.034
14	110.07–110.13	0.014–0.043	110.28–110.34	0.013–0.038	110.19–110.26	0.010–0.040
15	114.21–114.26	0.017–0.039	114.41–114.46	0.010–0.039	114.33–114.38	0.005–0.033
16	118.33–118.38	0.016–0.042	118.53–118.58	0.005–0.036	118.45–118.51	0.008–0.043
17	122.45–122.51	0.015–0.050	122.66–122.70	0.006–0.038	122.57–122.63	0.013–0.050
18	126.64–126.70	0.010–0.041	126.85–126.90	0.010–0.044	126.76–126.83	0.005–0.041
19	130.73–130.78	0.016–0.037	130.92–130.97	0.013–0.044	130.84–130.89	0.008–0.044
20	134.72–134.80	0.008–0.036	134.93–134.98	0.013–0.043	134.83–134.91	0.006–0.044
<b>vWA</b>						
11	151.38–151.45	0.015–0.052	151.76–151.80	0.010–0.045	151.65–151.76	0.019–0.048
12	155.39–155.45	0.021–0.048	155.78–155.83	0.005–0.044	155.69–155.77	0.010–0.049
13	159.46–159.52	0.018–0.049	159.85–159.90	0.010–0.037	159.76–159.83	0.008–0.045
14	163.68–163.74	0.024–0.055	164.07–164.12	0.010–0.041	163.97–164.05	0.010–0.053

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
15	167.62–167.69	0.026–0.047	168.01–168.06	0.010–0.045	167.93–168.00	0.010–0.050
16	171.66–171.73	0.021–0.049	172.05–172.11	0.010–0.050	171.98–172.06	0.010–0.059
17	175.71–175.77	0.017–0.053	176.10–176.15	0.005–0.046	176.05–176.10	0.013–0.047
18	179.72–179.77	0.014–0.054	180.10–180.18	0.005–0.034	180.03–180.11	0.013–0.060
19	183.76–183.83	0.029–0.054	184.16–184.22	0.008–0.046	184.12–184.17	0.000–0.057
20	187.79–187.87	0.028–0.054	188.20–188.26	0.005–0.056	188.12–188.21	0.015–0.057
21	191.80–191.86	0.023–0.048	192.20–192.26	0.005–0.040	192.13–192.21	0.013–0.067
22	195.83–195.88	0.021–0.055	196.23–196.29	0.006–0.051	196.14–196.24	0.013–0.051
23	199.77–199.84	0.027–0.048	200.20–200.25	0.010–0.046	200.12–200.19	0.010–0.062
24	204.14–204.20	0.027–0.058	204.54–204.60	0.013–0.049	204.48–204.53	0.013–0.057
<b>D16S539</b>						
5	226.98–227.05	0.029–0.055	227.54–227.60	0.008–0.043	227.44–227.50	0.029–0.077
8	239.10–239.18	0.037–0.067	239.67–239.73	0.013–0.043	239.59–239.64	0.034–0.079
9	243.17–243.24	0.029–0.070	243.73–243.79	0.000–0.046	243.60–243.70	0.025–0.090
10	247.21–247.27	0.041–0.078	247.77–247.83	0.013–0.050	247.67–247.75	0.017–0.068
11	251.26–251.33	0.025–0.066	251.84–251.87	0.010–0.045	251.72–251.80	0.026–0.081
12	255.30–255.36	0.043–0.061	255.88–255.93	0.015–0.045	255.76–255.83	0.029–0.075
13	259.33–259.39	0.040–0.074	259.92–259.98	0.008–0.053	259.81–259.88	0.017–0.088
14	263.39–263.45	0.038–0.077	263.96–264.02	0.008–0.057	263.86–263.94	0.021–0.077
15	267.42–267.48	0.037–0.077	268.01–268.07	0.005–0.056	267.89–267.99	0.025–0.094
<b>CSF1PO</b>						
6	281.95–282.03	0.035–0.077	282.56–282.63	0.010–0.057	282.49–282.58	0.029–0.107
7	285.99–286.07	0.042–0.076	286.58–286.65	0.016–0.045	286.53–286.61	0.024–0.099
8	290.02–290.10	0.047–0.068	290.61–290.69	0.016–0.047	290.57–290.65	0.021–0.090
9	294.04–294.13	0.042–0.071	294.65–294.72	0.008–0.042	294.60–294.69	0.027–0.104
10	298.07–298.15	0.039–0.069	298.68–298.75	0.010–0.050	298.65–298.73	0.028–0.087
11	302.10–302.18	0.033–0.074	302.72–302.79	0.006–0.059	302.66–302.75	0.040–0.081

**Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)**

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
12	306.12–306.20	0.032–0.075	306.77–306.83	0.013–0.061	306.72–306.79	0.029–0.093
13	310.15–310.23	0.042–0.069	310.78–310.85	0.013–0.054	310.76–310.83	0.024–0.095
14	314.17–314.27	0.045–0.081	314.83–314.89	0.013–0.051	314.78–314.87	0.033–0.091
15	318.19–318.28	0.021–0.072	318.85–318.93	0.010–0.065	318.84–318.91	0.033–0.085
<b>D6S1043</b>						
7	334.73–334.82	0.029–0.074	335.38–335.44	0.006–0.055	335.32–335.38	0.024–0.074
8	338.78–338.87	0.028–0.067	339.43–339.48	0.008–0.055	339.37–339.45	0.028–0.097
9	342.72–342.83	0.023–0.079	343.42–343.48	0.008–0.059	343.32–343.42	0.029–0.087
10	346.78–346.87	0.027–0.071	347.46–347.52	0.008–0.054	347.40–347.48	0.021–0.083
11	350.84–350.93	0.035–0.072	351.52–351.57	0.000–0.049	351.46–351.54	0.025–0.079
12	354.84–354.94	0.047–0.069	355.53–355.61	0.006–0.062	355.47–355.56	0.029–0.076
13	358.88–358.97	0.049–0.085	359.57–359.62	0.008–0.059	359.47–359.57	0.013–0.081
14	362.92–363.01	0.041–0.083	363.61–363.66	0.014–0.060	363.53–363.61	0.013–0.076
15	366.91–367.02	0.037–0.069	367.63–367.69	0.013–0.067	367.55–367.63	0.021–0.085
16	370.90–370.99	0.042–0.065	371.60–371.65	0.010–0.057	371.51–371.60	0.037–0.083
17	374.91–374.99	0.039–0.062	375.63–375.69	0.013–0.060	375.55–375.63	0.019–0.096
18	378.89–378.98	0.026–0.066	379.63–379.70	0.010–0.061	379.55–379.63	0.008–0.089
19	382.92–382.99	0.030–0.088	383.67–383.73	0.008–0.051	383.58–383.67	0.018–0.074
20	386.92–386.99	0.028–0.075	387.69–387.76	0.010–0.061	387.60–387.69	0.026–0.098
21	390.90–390.99	0.029–0.082	391.71–391.76	0.006–0.050	391.62–391.69	0.034–0.091
22	394.88–394.95	0.021–0.070	395.70–395.76	0.015–0.066	395.57–395.68	0.024–0.081
23	398.85–398.93	0.024–0.082	399.68–399.74	0.008–0.051	399.57–399.68	0.029–0.081
24	402.82–402.91	0.034–0.088	403.70–403.74	0.010–0.057	403.60–403.67	0.027–0.098
25	406.80–406.89	0.035–0.082	407.68–407.75	0.013–0.053	407.57–407.68	0.030–0.085
<b>Y Indel</b>						
1	81.20–81.26	0.019–0.047	81.43–81.47	0.005–0.051	81.33–81.38	0.008–0.036
2	86.50–86.56	0.023–0.041	86.73–86.78	0.008–0.045	86.62–86.70	0.010–0.042

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
<b>AMEL</b>						
X	102.45–102.50	0.014–0.044	102.53–102.57	0.008–0.034	102.44–102.50	0.013–0.041
Y	108.41–108.46	0.016–0.047	108.52–108.60	0.013–0.032	108.47–108.52	0.008–0.042
<b>D8S1179</b>						
5	114.53–114.58	0.019–0.035	114.64–114.68	0.008–0.038	114.57–114.62	0.005–0.054
6	118.63–118.69	0.014–0.052	118.74–118.79	0.005–0.039	118.67–118.74	0.006–0.044
7	122.74–122.79	0.009–0.040	122.84–122.89	0.006–0.046	122.75–122.83	0.010–0.047
8	126.84–126.89	0.009–0.039	126.94–126.99	0.010–0.047	126.88–126.93	0.005–0.045
9	130.93–131.00	0.008–0.045	131.03–131.08	0.010–0.040	130.97–131.05	0.010–0.042
10	135.02–135.08	0.014–0.041	135.12–135.16	0.010–0.047	135.04–135.11	0.005–0.061
11	139.11–139.16	0.012–0.047	139.19–139.24	0.008–0.058	139.13–139.18	0.000–0.048
12	143.20–143.27	0.015–0.039	143.29–143.34	0.013–0.040	143.23–143.29	0.013–0.047
13	147.39–147.44	0.012–0.035	147.47–147.51	0.010–0.039	147.40–147.47	0.008–0.052
14	151.47–151.53	0.016–0.046	151.52–151.58	0.006–0.039	151.48–151.54	0.010–0.054
15	155.54–155.59	0.010–0.037	155.58–155.64	0.005–0.039	155.54–155.60	0.010–0.048
16	159.61–159.65	0.016–0.032	159.65–159.71	0.010–0.048	159.58–159.65	0.010–0.046
17	163.66–163.72	0.010–0.032	163.70–163.75	0.010–0.041	163.66–163.71	0.006–0.059
18	167.73–167.78	0.015–0.035	167.76–167.82	0.013–0.047	167.72–167.76	0.010–0.045
19	171.79–171.86	0.008–0.027	171.82–171.86	0.010–0.044	171.77–171.83	0.010–0.051
<b>D21S11</b>						
24	182.85–182.91	0.033–0.079	183.23–183.28	0.006–0.049	183.17–183.25	0.005–0.070
24	184.88–184.94	0.028–0.061	185.26–185.31	0.008–0.056	185.23–185.28	0.013–0.071
25	186.88–186.95	0.037–0.068	187.28–187.32	0.005–0.056	187.23–187.28	0.019–0.062
26	190.94–190.98	0.034–0.076	191.32–191.37	0.008–0.044	191.27–191.33	0.017–0.072
27	194.98–195.04	0.042–0.065	195.38–195.42	0.008–0.047	195.31–195.40	0.014–0.077
28	198.99–199.05	0.035–0.077	199.37–199.43	0.010–0.044	199.34–199.40	0.013–0.067
28	200.99–201.03	0.039–0.061	201.38–201.43	0.010–0.045	201.34–201.39	0.008–0.074



**Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)**

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
29	203.00–203.05	0.045–0.077	203.39–203.44	0.008–0.052	203.37–203.42	0.005–0.067
29	205.06–205.10	0.042–0.079	205.44–205.50	0.015–0.047	205.40–205.47	0.021–0.078
30	207.05–207.10	0.039–0.074	207.45–207.51	0.013–0.045	207.39–207.47	0.017–0.082
30	209.05–209.11	0.042–0.072	209.45–209.49	0.010–0.043	209.39–209.47	0.021–0.067
31	211.08–211.14	0.032–0.067	211.49–211.55	0.008–0.045	211.43–211.51	0.010–0.072
31	213.08–213.15	0.039–0.073	213.49–213.52	0.010–0.043	213.44–213.50	0.014–0.081
32	215.12–215.17	0.042–0.074	215.52–215.57	0.005–0.040	215.47–215.54	0.010–0.069
32	217.10–217.16	0.050–0.069	217.52–217.56	0.005–0.048	217.44–217.54	0.010–0.069
33	219.14–219.22	0.048–0.075	219.55–219.61	0.006–0.039	219.52–219.59	0.017–0.080
33	221.10–221.16	0.044–0.071	221.50–221.57	0.010–0.050	221.48–221.54	0.020–0.085
34	223.23–223.29	0.035–0.078	223.65–223.69	0.013–0.040	223.59–223.66	0.022–0.075
34	225.17–225.23	0.043–0.074	225.58–225.64	0.010–0.042	225.54–225.61	0.021–0.090
35	227.26–227.30	0.030–0.080	227.66–227.71	0.005–0.046	227.62–227.68	0.024–0.077
35	229.20–229.25	0.042–0.073	229.61–229.67	0.005–0.048	229.57–229.64	0.022–0.076
36	231.20–231.26	0.037–0.082	231.63–231.69	0.010–0.051	231.58–231.65	0.017–0.075
37	235.28–235.34	0.042–0.076	235.71–235.75	0.006–0.051	235.65–235.73	0.030–0.070
38	239.25–239.32	0.040–0.082	239.67–239.75	0.008–0.045	239.62–239.71	0.028–0.072
<b>D18S51</b>						
7	261.02–261.08	0.037–0.075	261.35–261.42	0.005–0.044	261.32–261.38	0.037–0.064
9	269.17–269.23	0.028–0.068	269.49–269.56	0.010–0.051	269.44–269.51	0.010–0.068
10	273.24–273.30	0.034–0.062	273.56–273.62	0.013–0.045	273.52–273.59	0.013–0.071
10	275.24–275.30	0.034–0.079	275.56–275.61	0.010–0.041	275.51–275.60	0.024–0.060
11	277.31–277.37	0.025–0.063	277.63–277.69	0.010–0.046	277.58–277.65	0.015–0.062
12	281.38–281.45	0.031–0.073	281.70–281.75	0.008–0.056	281.64–281.72	0.015–0.078
13	285.46–285.53	0.037–0.069	285.77–285.81	0.010–0.051	285.69–285.79	0.016–0.063
13	287.48–287.54	0.027–0.075	287.77–287.82	0.017–0.051	287.71–287.79	0.010–0.062
14	289.53–289.60	0.034–0.058	289.82–289.89	0.013–0.044	289.80–289.87	0.013–0.068

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
14	291.56–291.61	0.032–0.069	291.82–291.89	0.013–0.051	291.79–291.85	0.005–0.070
15	293.62–293.68	0.020–0.057	293.91–293.96	0.008–0.046	293.85–293.93	0.010–0.064
16	297.70–297.77	0.035–0.064	297.97–298.03	0.010–0.059	297.92–298.00	0.010–0.059
17	301.76–301.83	0.029–0.071	302.03–302.09	0.017–0.051	302.01–302.08	0.019–0.062
18	305.85–305.90	0.025–0.078	306.11–306.18	0.010–0.055	306.08–306.14	0.021–0.053
19	309.94–309.98	0.029–0.061	310.16–310.22	0.016–0.051	310.15–310.21	0.008–0.067
20	314.00–314.06	0.021–0.054	314.25–314.29	0.013–0.045	314.21–314.28	0.010–0.059
21	318.10–318.13	0.021–0.057	318.31–318.37	0.010–0.043	318.26–318.37	0.008–0.054
22	322.22–322.28	0.015–0.055	322.44–322.50	0.005–0.053	322.42–322.49	0.010–0.062
23	326.22–326.28	0.025–0.064	326.43–326.51	0.013–0.057	326.42–326.49	0.015–0.076
24	330.27–330.34	0.022–0.067	330.51–330.56	0.010–0.041	330.49–330.57	0.013–0.073
25	334.34–334.41	0.018–0.062	334.56–334.63	0.008–0.056	334.56–334.63	0.010–0.059
26	338.44–338.49	0.022–0.055	338.63–338.69	0.005–0.050	338.63–338.72	0.010–0.055
27	342.48–342.56	0.024–0.068	342.71–342.76	0.006–0.052	342.72–342.79	0.014–0.059
<b>D5S818</b>						
7	370.14–370.21	0.037–0.079	370.66–370.71	0.015–0.068	370.60–370.70	0.006–0.077
8	374.15–374.21	0.029–0.059	374.68–374.75	0.010–0.064	374.65–374.73	0.013–0.079
9	378.14–378.20	0.019–0.063	378.71–378.76	0.010–0.074	378.67–378.75	0.018–0.076
10	382.14–382.20	0.016–0.068	382.71–382.79	0.010–0.065	382.69–382.77	0.013–0.076
11	386.13–386.20	0.024–0.072	386.74–386.80	0.008–0.057	386.71–386.78	0.024–0.090
12	390.12–390.19	0.037–0.072	390.75–390.82	0.013–0.061	390.73–390.80	0.021–0.082
13	394.10–394.18	0.038–0.102	394.76–394.83	0.015–0.067	394.72–394.82	0.032–0.087
14	398.10–398.16	0.032–0.071	398.77–398.82	0.010–0.068	398.74–398.82	0.031–0.091
15	402.08–402.14	0.026–0.073	402.77–402.84	0.012–0.058	402.73–402.81	0.017–0.095
16	406.06–406.12	0.032–0.110	406.78–406.85	0.013–0.075	406.73–406.81	0.017–0.085
17	410.04–410.10	0.044–0.114	410.78–410.85	0.008–0.079	410.72–410.82	0.021–0.100
18	414.00–414.08	0.043–0.103	414.77–414.85	0.017–0.057	414.73–414.80	0.037–0.106

**Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)**

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
<b>D2S441</b>						
8	83.24–83.27	0.008–0.034	83.25–83.30	0.005–0.050	83.19–83.25	0.005–0.036
9	87.34–87.37	0.012–0.031	87.37–87.42	0.006–0.038	87.31–87.40	0.008–0.036
10	91.40–91.42	0.010–0.034	91.45–91.50	0.005–0.041	91.40–91.45	0.008–0.035
11	95.52–95.56	0.014–0.046	95.60–95.65	0.006–0.046	95.56–95.61	0.013–0.038
11	98.49–98.53	0.015–0.032	98.57–98.63	0.008–0.039	98.54–98.58	0.008–0.048
12	99.59–99.64	0.015–0.037	99.69–99.73	0.008–0.041	99.65–99.70	0.008–0.035
13	103.52–103.55	0.018–0.036	103.62–103.68	0.013–0.036	103.58–103.64	0.013–0.045
14	107.59–107.63	0.015–0.049	107.71–107.76	0.014–0.038	107.67–107.71	0.008–0.046
15	111.64–111.69	0.015–0.045	111.78–111.82	0.014–0.045	111.75–111.80	0.006–0.039
16	115.73–115.78	0.010–0.033	115.85–115.90	0.010–0.034	115.81–115.89	0.010–0.044
17	119.88–119.93	0.014–0.033	120.02–120.07	0.013–0.045	120.01–120.05	0.013–0.039
<b>D19S433</b>						
5	141.04–141.09	0.016–0.043	141.42–141.46	0.006–0.050	141.37–141.44	0.021–0.059
6	145.02–145.07	0.019–0.059	145.41–145.45	0.008–0.038	145.39–145.45	0.013–0.071
7	148.99–149.05	0.015–0.048	149.39–149.44	0.014–0.051	149.38–149.44	0.010–0.063
8	152.95–153.02	0.014–0.056	153.38–153.42	0.010–0.046	153.33–153.42	0.018–0.069
9	156.91–156.98	0.026–0.055	157.35–157.41	0.008–0.042	157.35–157.40	0.014–0.073
10	160.86–160.94	0.029–0.051	161.33–161.38	0.010–0.041	161.32–161.38	0.013–0.065
11	164.82–164.89	0.017–0.048	165.30–165.36	0.010–0.052	165.28–165.36	0.017–0.070
12	168.78–168.85	0.029–0.054	169.26–169.32	0.008–0.052	169.24–169.33	0.024–0.084
12	170.89–170.97	0.032–0.059	171.38–171.44	0.006–0.046	171.37–171.45	0.005–0.076
13	172.74–172.81	0.018–0.060	173.24–173.29	0.005–0.048	173.22–173.30	0.027–0.080
13	174.72–174.80	0.025–0.062	175.22–175.28	0.010–0.045	175.21–175.30	0.027–0.079
14	176.68–176.76	0.030–0.065	177.20–177.26	0.008–0.054	177.16–177.27	0.028–0.082
14	178.65–178.74	0.026–0.069	179.17–179.25	0.005–0.037	179.16–179.27	0.018–0.088
15	180.62–180.70	0.029–0.070	181.14–181.21	0.005–0.039	181.11–181.23	0.024–0.098

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
15	182.63–182.69	0.028–0.064	183.15–183.22	0.006–0.046	183.15–183.22	0.028–0.088
16	184.56–184.66	0.034–0.060	185.10–185.17	0.006–0.046	185.07–185.19	0.031–0.077
16	186.57–186.66	0.028–0.063	187.11–187.17	0.005–0.043	187.13–187.20	0.029–0.083
17	188.49–188.59	0.028–0.070	189.06–189.13	0.005–0.053	189.05–189.15	0.032–0.087
17	190.53–190.62	0.031–0.064	191.09–191.15	0.005–0.062	191.07–191.16	0.030–0.096
18	194.56–194.64	0.032–0.060	195.12–195.19	0.008–0.032	195.12–195.20	0.037–0.077
19	198.49–198.58	0.031–0.062	199.08–199.14	0.010–0.051	199.08–199.17	0.028–0.105
<b>FGA</b>						
13	207.89–207.92	0.012–0.039	207.88–207.93	0.008–0.052	207.83–207.89	0.010–0.042
14	211.94–211.97	0.013–0.030	211.93–211.97	0.010–0.051	211.88–211.94	0.013–0.055
15	215.99–216.02	0.011–0.035	215.96–216.02	0.008–0.051	215.93–215.97	0.005–0.049
16	220.05–220.09	0.016–0.032	220.01–220.07	0.008–0.047	219.97–220.04	0.010–0.046
17	224.11–224.15	0.015–0.039	224.05–224.11	0.010–0.043	224.01–224.08	0.010–0.040
18	228.16–228.20	0.013–0.031	228.11–228.17	0.006–0.053	228.07–228.12	0.005–0.049
19	232.22–232.27	0.013–0.033	232.15–232.20	0.010–0.042	232.12–232.17	0.010–0.048
20	236.27–236.33	0.010–0.041	236.20–236.25	0.010–0.046	236.17–236.22	0.005–0.043
21	240.33–240.38	0.008–0.052	240.25–240.30	0.010–0.045	240.21–240.27	0.006–0.039
22	244.39–244.44	0.012–0.034	244.30–244.35	0.000–0.045	244.25–244.33	0.012–0.041
23	248.46–248.50	0.012–0.044	248.34–248.40	0.013–0.038	248.31–248.36	0.008–0.050
24	252.51–252.56	0.014–0.040	252.39–252.45	0.010–0.046	252.38–252.43	0.010–0.044
25	256.58–256.62	0.016–0.033	256.46–256.51	0.014–0.044	256.41–256.49	0.010–0.042
26	260.63–260.68	0.014–0.046	260.48–260.54	0.005–0.052	260.46–260.52	0.010–0.050
26	262.66–262.71	0.014–0.063	262.53–262.59	0.010–0.050	262.49–262.55	0.005–0.041
27	264.66–264.70	0.012–0.040	264.52–264.56	0.010–0.041	264.48–264.53	0.010–0.048
28	268.72–268.76	0.010–0.046	268.58–268.61	0.010–0.048	268.53–268.59	0.008–0.045
29	272.78–272.83	0.017–0.063	272.63–272.69	0.013–0.038	272.61–272.66	0.005–0.049
30	276.86–276.92	0.011–0.048	276.71–276.78	0.008–0.045	276.68–276.73	0.006–0.045

**Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)**

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
30	278.69–278.75	0.012–0.044	278.55–278.60	0.010–0.053	278.51–278.58	0.006–0.048
31	282.76–282.81	0.012–0.034	282.59–282.67	0.008–0.045	282.55–282.63	0.008–0.048
32	286.83–286.88	0.018–0.038	286.68–286.73	0.010–0.042	286.62–286.69	0.013–0.051
33	290.89–290.94	0.014–0.036	290.71–290.77	0.013–0.041	290.67–290.75	0.008–0.049
42	327.62–327.67	0.014–0.050	327.39–327.43	0.010–0.053	327.38–327.46	0.008–0.057
43	331.67–331.73	0.006–0.061	331.44–331.49	0.010–0.049	331.42–331.51	0.015–0.063
44	335.72–335.79	0.015–0.046	335.50–335.57	0.010–0.046	335.53–335.59	0.013–0.062
45	339.80–339.86	0.019–0.048	339.57–339.63	0.000–0.050	339.59–339.65	0.010–0.068
46	343.77–343.84	0.013–0.054	343.52–343.59	0.005–0.057	343.54–343.61	0.005–0.077
47	347.84–347.90	0.013–0.047	347.57–347.64	0.013–0.041	347.59–347.67	0.013–0.067
48	351.92–351.98	0.012–0.043	351.66–351.72	0.010–0.048	351.67–351.75	0.006–0.076
49	355.98–356.04	0.014–0.044	355.71–355.77	0.000–0.045	355.74–355.80	0.012–0.060
50	360.03–360.07	0.018–0.054	359.73–359.80	0.013–0.050	359.75–359.82	0.017–0.079
51	364.08–364.14	0.018–0.078	363.80–363.85	0.014–0.056	363.82–363.88	0.017–0.073
<b>D10S1248</b>						
8	371.77–371.82	0.023–0.063	372.17–372.23	0.010–0.057	372.12–372.21	0.010–0.071
9	375.73–375.80	0.023–0.061	376.16–376.23	0.010–0.059	376.13–376.20	0.017–0.069
10	379.73–379.79	0.031–0.058	380.16–380.23	0.010–0.047	380.14–380.19	0.016–0.056
11	383.68–383.76	0.024–0.060	384.12–384.19	0.008–0.054	384.09–384.16	0.010–0.068
12	387.70–387.76	0.020–0.056	388.15–388.22	0.012–0.054	388.13–388.18	0.021–0.058
13	391.69–391.75	0.019–0.163	392.14–392.20	0.013–0.062	392.11–392.17	0.010–0.072
14	395.65–395.72	0.028–0.088	396.10–396.17	0.013–0.054	396.07–396.15	0.008–0.076
15	399.64–399.72	0.028–0.128	400.11–400.17	0.013–0.054	400.07–400.14	0.017–0.074
16	403.61–403.67	0.024–0.062	404.10–404.19	0.013–0.058	404.07–404.14	0.017–0.077
17	407.63–407.68	0.025–0.081	408.13–408.18	0.005–0.057	408.08–408.14	0.017–0.085
18	411.56–411.60	0.019–0.115	412.04–412.12	0.008–0.052	412.00–412.09	0.014–0.067
19	415.49–415.57	0.021–0.102	416.02–416.10	0.015–0.056	416.00–416.05	0.017–0.080

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
<b>D22S1045</b>						
8	88.43–88.47	0.012–0.037	88.46–88.53	0.005–0.039	88.40–88.45	0.005–0.052
9	91.42–91.46	0.013–0.037	91.46–91.54	0.005–0.042	91.41–91.46	0.010–0.042
10	94.42–94.46	0.010–0.030	94.50–94.54	0.005–0.042	94.43–94.47	0.012–0.039
11	97.43–97.46	0.013–0.038	97.49–97.56	0.008–0.045	97.43–97.49	0.005–0.054
12	100.42–100.45	0.010–0.042	100.51–100.56	0.006–0.043	100.45–100.50	0.005–0.036
13	103.41–103.46	0.012–0.038	103.50–103.57	0.008–0.040	103.44–103.49	0.013–0.041
14	106.40–106.44	0.008–0.039	106.51–106.55	0.015–0.045	106.43–106.48	0.014–0.041
15	109.39–109.44	0.013–0.044	109.51–109.57	0.013–0.039	109.44–109.50	0.010–0.043
16	112.38–112.42	0.017–0.038	112.50–112.54	0.010–0.039	112.43–112.49	0.000–0.037
17	115.37–115.40	0.008–0.028	115.49–115.56	0.010–0.044	115.41–115.49	0.013–0.040
18	118.35–118.41	0.012–0.032	118.50–118.55	0.013–0.053	118.41–118.48	0.008–0.041
19	121.34–121.40	0.010–0.040	121.48–121.53	0.005–0.042	121.41–121.46	0.006–0.042
<b>D1S1656</b>						
9	146.83–146.86	0.011–0.039	146.87–146.93	0.010–0.043	146.86–146.91	0.005–0.041
10	150.72–150.76	0.015–0.034	150.77–150.83	0.010–0.039	150.78–150.81	0.017–0.043
11	154.72–154.77	0.010–0.038	154.78–154.83	0.008–0.044	154.77–154.84	0.014–0.050
12	158.81–158.85	0.014–0.031	158.86–158.92	0.010–0.042	158.88–158.91	0.013–0.048
13	162.74–162.76	0.017–0.034	162.80–162.85	0.014–0.037	162.80–162.85	0.010–0.043
14	166.81–166.85	0.012–0.038	166.89–166.93	0.013–0.048	166.87–166.94	0.010–0.051
14	169.77–169.80	0.010–0.034	169.85–169.89	0.008–0.043	169.85–169.89	0.010–0.047
15	170.79–170.83	0.010–0.047	170.87–170.92	0.006–0.036	170.86–170.93	0.000–0.051
15	173.75–173.79	0.014–0.035	173.84–173.90	0.005–0.042	173.83–173.89	0.006–0.043
16	174.80–174.85	0.014–0.035	174.88–174.93	0.000–0.036	174.88–174.94	0.008–0.053
16	177.75–177.80	0.016–0.040	177.83–177.89	0.008–0.044	177.84–177.89	0.010–0.048
17	178.88–178.92	0.012–0.064	178.95–179.00	0.008–0.040	178.95–179.00	0.013–0.053
17	181.75–181.79	0.009–0.043	181.84–181.89	0.005–0.039	181.82–181.89	0.014–0.044

**Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)**

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
18	182.78–182.82	0.010–0.034	182.87–182.92	0.005–0.050	182.85–182.92	0.005–0.051
18	185.75–185.78	0.012–0.042	185.84–185.89	0.005–0.041	185.85–185.90	0.010–0.040
19	189.77–189.80	0.017–0.036	189.88–189.93	0.006–0.043	189.87–189.93	0.006–0.052
20	193.82–193.86	0.014–0.035	193.93–193.98	0.005–0.033	193.92–193.98	0.008–0.059
<b>D13S317</b>						
5	201.31–201.35	0.020–0.036	201.46–201.52	0.010–0.047	201.45–201.50	0.013–0.053
6	205.33–205.37	0.015–0.031	205.47–205.52	0.015–0.038	205.46–205.51	0.005–0.045
7	209.35–209.39	0.011–0.039	209.49–209.54	0.010–0.050	209.48–209.53	0.017–0.056
8	213.37–213.42	0.015–0.033	213.49–213.55	0.005–0.039	213.50–213.55	0.014–0.049
9	217.42–217.46	0.014–0.036	217.53–217.58	0.005–0.039	217.50–217.56	0.005–0.048
10	221.43–221.48	0.019–0.036	221.55–221.60	0.010–0.054	221.53–221.60	0.006–0.045
11	225.48–225.50	0.018–0.038	225.58–225.63	0.010–0.043	225.57–225.62	0.017–0.048
12	229.49–229.54	0.014–0.039	229.60–229.66	0.008–0.049	229.59–229.64	0.010–0.043
13	233.54–233.58	0.012–0.044	233.62–233.68	0.008–0.043	233.63–233.66	0.015–0.051
14	237.55–237.61	0.012–0.034	237.66–237.72	0.005–0.044	237.64–237.69	0.006–0.054
15	241.59–241.64	0.005–0.041	241.69–241.74	0.006–0.048	241.67–241.72	0.010–0.053
16	245.64–245.70	0.015–0.043	245.72–245.78	0.008–0.042	245.71–245.76	0.015–0.066
<b>D7S820</b>						
6	259.89–259.94	0.010–0.048	260.03–260.07	0.008–0.049	259.99–260.05	0.008–0.046
7	263.90–263.96	0.008–0.045	264.06–264.11	0.013–0.042	264.00–264.07	0.014–0.049
8	267.92–267.97	0.018–0.044	268.08–268.13	0.010–0.048	268.05–268.09	0.010–0.056
9	271.95–271.99	0.009–0.045	272.10–272.15	0.017–0.045	272.07–272.12	0.010–0.050
10	275.96–276.01	0.017–0.042	276.13–276.19	0.010–0.047	276.10–276.16	0.010–0.061
11	279.97–280.02	0.012–0.057	280.16–280.22	0.010–0.047	280.10–280.19	0.014–0.052
12	284.00–284.04	0.017–0.042	284.20–284.25	0.012–0.039	284.14–284.21	0.012–0.056
13	288.02–288.08	0.012–0.040	288.21–288.28	0.014–0.042	288.18–288.24	0.008–0.061
14	292.05–292.09	0.013–0.045	292.24–292.30	0.013–0.048	292.21–292.27	0.010–0.062

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
15	296.07–296.11	0.017–0.040	296.26–296.33	0.010–0.052	296.24–296.28	0.010–0.053
<b>Penta E</b>						
5	310.56–310.66	0.075–0.110	311.23–311.28	0.010–0.059	311.20–311.26	0.044–0.096
6	315.69–315.78	0.068–0.113	316.34–316.39	0.014–0.055	316.30–316.38	0.014–0.097
7	320.72–320.82	0.073–0.121	321.41–321.47	0.005–0.051	321.39–321.46	0.028–0.099
8	325.78–325.88	0.060–0.108	326.44–326.50	0.010–0.057	326.42–326.48	0.029–0.096
9	330.82–330.91	0.060–0.109	331.48–331.54	0.008–0.052	331.46–331.54	0.025–0.090
10	335.88–335.97	0.071–0.103	336.53–336.59	0.010–0.058	336.53–336.60	0.040–0.097
11	340.87–341.01	0.064–0.117	341.58–341.66	0.005–0.056	341.59–341.65	0.040–0.094
12	345.96–346.05	0.078–0.118	346.63–346.69	0.013–0.070	346.64–346.71	0.036–0.081
13	351.02–351.12	0.065–0.120	351.69–351.75	0.005–0.055	351.70–351.78	0.017–0.091
14	356.06–356.18	0.075–0.123	356.74–356.80	0.018–0.066	356.74–356.83	0.030–0.090
15	361.13–361.23	0.093–0.137	361.80–361.86	0.013–0.055	361.78–361.87	0.033–0.090
16	366.21–366.29	0.073–0.122	366.83–366.91	0.013–0.067	366.82–366.91	0.019–0.082
17	371.24–371.35	0.069–0.117	371.90–371.98	0.013–0.070	371.87–371.96	0.021–0.080
18	376.29–376.36	0.066–0.116	376.96–377.02	0.019–0.068	376.93–377.00	0.030–0.083
19	381.31–381.41	0.070–0.104	382.00–382.07	0.014–0.059	381.98–382.03	0.012–0.083
20	386.35–386.43	0.069–0.108	387.04–387.09	0.010–0.070	387.01–387.10	0.031–0.109
21	391.38–391.46	0.068–0.106	392.07–392.15	0.008–0.056	392.05–392.14	0.005–0.087
22	396.39–396.48	0.064–0.108	397.11–397.17	0.012–0.067	397.07–397.16	0.017–0.095
23	401.41–401.51	0.071–0.111	402.15–402.21	0.010–0.054	402.11–402.21	0.030–0.101
24	406.40–406.52	0.077–0.120	407.17–407.23	0.005–0.052	407.14–407.22	0.013–0.092
25	411.44–411.54	0.081–0.109	412.20–412.25	0.008–0.051	412.16–412.24	0.029–0.083
26	416.45–416.52	0.076–0.113	417.22–417.27	0.017–0.067	417.21–417.28	0.026–0.095
<b>Penta D</b>						
2	78.64–78.68	0.008–0.055	78.99–79.05	0.010–0.048	78.87–78.95	0.010–0.061
3	83.57–83.62	0.013–0.057	83.98–84.05	0.005–0.049	83.86–83.94	0.017–0.063



**Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)**

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
5	91.30–91.35	0.023–0.053	91.78–91.86	0.010–0.037	91.65–91.74	0.018–0.062
6	96.21–96.27	0.023–0.067	96.71–96.79	0.010–0.041	96.62–96.69	0.022–0.064
7	101.32–101.37	0.023–0.060	101.85–101.92	0.008–0.035	101.72–101.81	0.013–0.059
8	105.99–106.05	0.028–0.060	106.60–106.65	0.012–0.044	106.47–106.57	0.022–0.077
9	110.88–110.96	0.028–0.063	111.52–111.57	0.010–0.033	111.38–111.48	0.025–0.071
10	115.91–115.96	0.025–0.060	116.55–116.60	0.005–0.041	116.44–116.51	0.021–0.073
11	120.62–120.69	0.038–0.075	121.30–121.38	0.008–0.041	121.17–121.28	0.029–0.076
12	125.49–125.57	0.035–0.073	126.21–126.27	0.017–0.042	126.07–126.19	0.030–0.077
13	130.45–130.53	0.031–0.078	131.21–131.28	0.010–0.038	131.06–131.18	0.031–0.081
14	135.19–135.27	0.033–0.077	135.95–136.04	0.008–0.049	135.82–135.94	0.029–0.094
15	140.05–140.13	0.038–0.073	140.85–140.93	0.010–0.043	140.73–140.85	0.039–0.092
16	145.02–145.10	0.046–0.077	145.85–145.90	0.008–0.042	145.70–145.81	0.029–0.098
17	149.75–149.84	0.048–0.081	150.61–150.67	0.010–0.061	150.44–150.59	0.043–0.102
<b>TH01</b>						
4	169.86–169.92	0.029–0.060	170.35–170.40	0.005–0.045	170.30–170.38	0.017–0.069
5	173.87–173.94	0.018–0.055	174.34–174.42	0.000–0.047	174.30–174.39	0.013–0.068
6	177.86–177.93	0.023–0.060	178.36–178.41	0.008–0.048	178.31–178.38	0.017–0.076
7	181.86–181.94	0.030–0.053	182.36–182.40	0.005–0.036	182.32–182.40	0.021–0.075
8	185.86–185.94	0.022–0.060	186.35–186.43	0.005–0.043	186.33–186.40	0.016–0.078
9	189.87–189.95	0.035–0.059	190.39–190.43	0.008–0.051	190.34–190.41	0.024–0.076
9	192.94–193.01	0.028–0.069	193.43–193.49	0.010–0.050	193.41–193.48	0.019–0.080
10	193.88–193.95	0.031–0.062	194.37–194.44	0.005–0.051	194.35–194.42	0.015–0.074
11	197.87–197.96	0.027–0.060	198.40–198.47	0.006–0.046	198.37–198.44	0.021–0.078
12	201.89–201.97	0.029–0.070	202.39–202.48	0.013–0.050	202.37–202.44	0.034–0.071
13	208.95–209.03	0.036–0.065	209.45–209.52	0.010–0.048	209.43–209.52	0.025–0.081
<b>D12S391</b>						
14	218.79–218.84	0.014–0.034	218.81–218.88	0.006–0.034	218.81–218.87	0.010–0.047

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
15	222.81–222.84	0.016–0.041	222.81–222.87	0.010–0.046	222.82–222.87	0.012–0.042
16	226.76–226.80	0.010–0.033	226.76–226.83	0.012–0.050	226.77–226.82	0.015–0.056
17	230.84–230.89	0.013–0.036	230.84–230.90	0.008–0.054	230.85–230.90	0.008–0.064
18	234.88–234.92	0.015–0.036	234.87–234.92	0.005–0.058	234.87–234.91	0.010–0.043
19	238.94–238.99	0.015–0.037	238.94–238.99	0.010–0.047	238.92–238.98	0.005–0.048
19	241.95–242.00	0.010–0.076	241.93–241.98	0.005–0.043	241.93–241.97	0.010–0.055
20	242.89–242.95	0.012–0.043	242.91–242.95	0.010–0.052	242.89–242.95	0.008–0.044
21	247.01–247.06	0.010–0.043	246.99–247.04	0.010–0.041	246.97–247.04	0.005–0.046
22	251.12–251.17	0.012–0.038	251.09–251.15	0.014–0.041	251.08–251.13	0.013–0.050
23	255.12–255.18	0.011–0.033	255.10–255.15	0.010–0.040	255.09–255.15	0.013–0.042
24	259.13–259.19	0.018–0.034	259.11–259.16	0.010–0.042	259.09–259.15	0.010–0.033
25	263.16–263.22	0.017–0.035	263.13–263.19	0.013–0.053	263.12–263.18	0.008–0.046
26	267.23–267.27	0.015–0.048	267.19–267.23	0.006–0.038	267.16–267.22	0.005–0.043
27	271.22–271.27	0.012–0.041	271.18–271.24	0.013–0.046	271.16–271.23	0.010–0.045
<b>D2S1338</b>						
11	280.86–280.95	0.042–0.072	281.45–281.48	0.008–0.058	281.40–281.48	0.040–0.085
12	284.85–284.94	0.031–0.069	285.45–285.50	0.017–0.043	285.41–285.48	0.041–0.093
13	288.87–288.94	0.031–0.075	289.42–289.51	0.013–0.057	289.41–289.48	0.033–0.089
14	292.86–292.96	0.024–0.065	293.43–293.48	0.010–0.062	293.40–293.50	0.021–0.088
15	296.82–296.91	0.025–0.061	297.35–297.39	0.006–0.048	297.31–297.39	0.021–0.089
16	300.87–300.94	0.034–0.067	301.39–301.45	0.015–0.053	301.37–301.44	0.029–0.086
17	304.87–304.94	0.033–0.063	305.40–305.46	0.010–0.054	305.39–305.46	0.021–0.088
18	308.91–308.98	0.029–0.070	309.41–309.48	0.014–0.053	309.41–309.47	0.029–0.082
19	312.91–312.98	0.034–0.058	313.44–313.50	0.013–0.053	313.41–313.48	0.039–0.091
20	316.92–317.00	0.020–0.058	317.45–317.51	0.013–0.060	317.43–317.49	0.021–0.088
21	320.97–321.05	0.021–0.063	321.48–321.54	0.005–0.051	321.47–321.54	0.027–0.084
22	324.97–325.06	0.033–0.070	325.48–325.53	0.010–0.053	325.46–325.53	0.015–0.070

Table 7 Precision results of multiple runs of the Verifiler Plus Allelic Ladder (continued)

Allele	3130xI		3500		3500xL	
	Mean	Standard deviation	Mean	Standard deviation	Mean	Standard deviation
23	328.98–329.06	0.024–0.060	329.49–329.55	0.013–0.073	329.48–329.58	0.015–0.088
24	333.00–333.08	0.024–0.061	333.53–333.58	0.010–0.054	333.51–333.61	0.015–0.075
25	337.01–337.10	0.028–0.063	337.53–337.58	0.010–0.066	337.54–337.63	0.015–0.080
26	341.06–341.13	0.026–0.063	341.54–341.60	0.005–0.059	341.56–341.64	0.015–0.078
27	345.14–345.22	0.028–0.067	345.61–345.68	0.008–0.039	345.61–345.69	0.006–0.079
28	349.38–349.46	0.021–0.062	349.77–349.81	0.006–0.055	349.80–349.86	0.010–0.074
<b>TPOX</b>						
5	364.37–364.50	0.073–0.107	365.42–365.42	0.013–0.013	365.34–365.44	0.043–0.119
6	368.42–368.52	0.061–0.110	369.45–369.45	0.010–0.010	369.37–369.50	0.040–0.112
7	372.46–372.59	0.054–0.102	373.51–373.51	0.008–0.008	373.45–373.54	0.040–0.133
8	376.44–376.57	0.053–0.087	377.49–377.49	0.010–0.010	377.44–377.54	0.044–0.110
9	380.42–380.54	0.043–0.090	381.52–381.52	0.014–0.014	381.43–381.55	0.050–0.113
10	384.44–384.56	0.037–0.093	385.53–385.53	0.005–0.005	385.47–385.58	0.046–0.142
11	388.43–388.55	0.054–0.092	389.57–389.57	0.010–0.010	389.51–389.58	0.055–0.128
12	392.41–392.53	0.046–0.098	393.59–393.59	0.015–0.015	393.50–393.60	0.065–0.121
13	396.42–396.53	0.042–0.095	397.59–397.59	0.010–0.010	397.52–397.63	0.066–0.124
14	400.41–400.53	0.050–0.102	401.61–401.61	0.013–0.013	401.50–401.66	0.061–0.127
15	404.39–404.51	0.057–0.098	405.63–405.63	0.008–0.008	405.51–405.64	0.047–0.141

## 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.

### 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 observation

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.
- Smaller alleles typically show a lower level of stutter relative to the longer alleles in each locus.
- Each allele in a locus displays a consistent stutter percentage.
- Peaks in the stutter position that are above the stutter filter percentage specified in the software are not filtered. Peaks in the stutter position that have not been filtered and remain labeled can be further evaluated.
- 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.
- The magnitude and/or variability of stutter may increase with low DNA input amounts.

The stutter observed in the population study that are one repeat unit away from the alleles recorded are shown in Figure 23 through Figure 27. All data were generated on the 3500xL Genetic Analyzer.

The stutter filter settings that are derived from these data are listed in “Stutter filter settings provided with the GeneMapper™ ID-X Software” on page 81.

Plus-stutter was regularly observed and was more significant in trinucleotide repeat-containing loci (see “Plus-stutter peaks at the D22S1045 locus” on page 80).

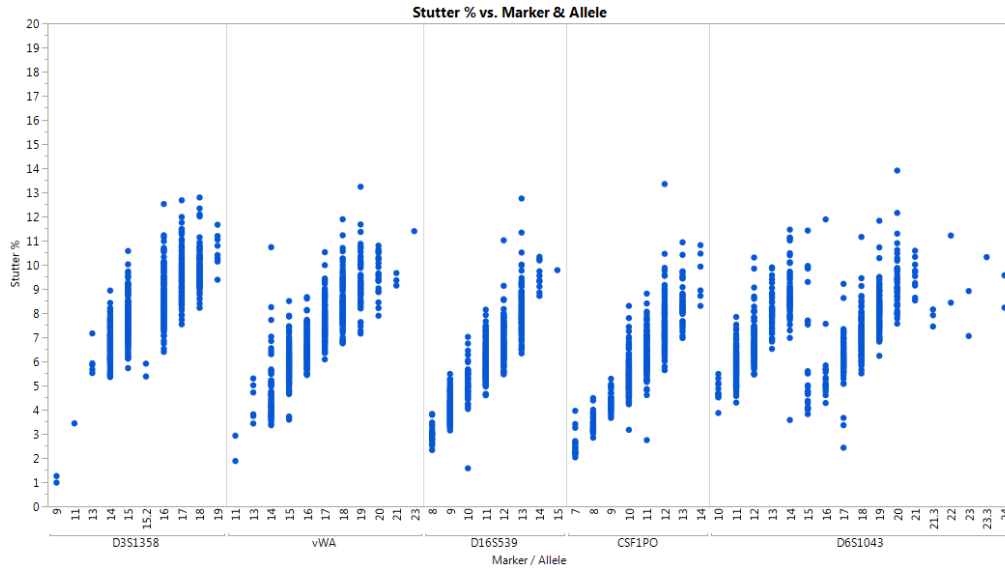


Figure 23 Stutter percentages for the Verifiler™ Plus kit FAM™ dye (blue) channel loci: D3S1358, vWA, D16S539, CSF1PO, and D6S1043.

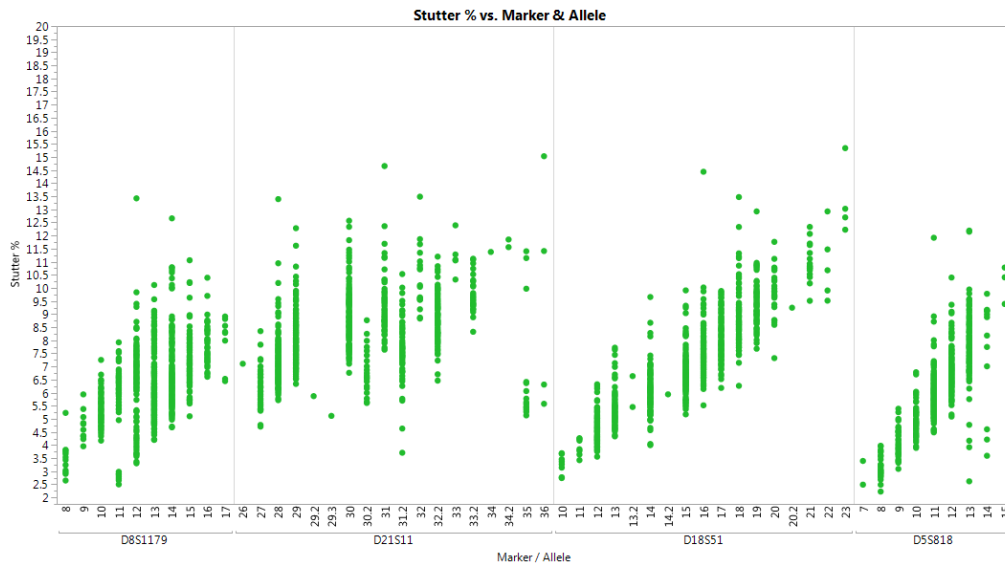


Figure 24 Stutter percentages for the Verifiler™ Plus kit VIC™ dye (green) channel loci: D8S1179, D21S11, D18S51, and D5S818.

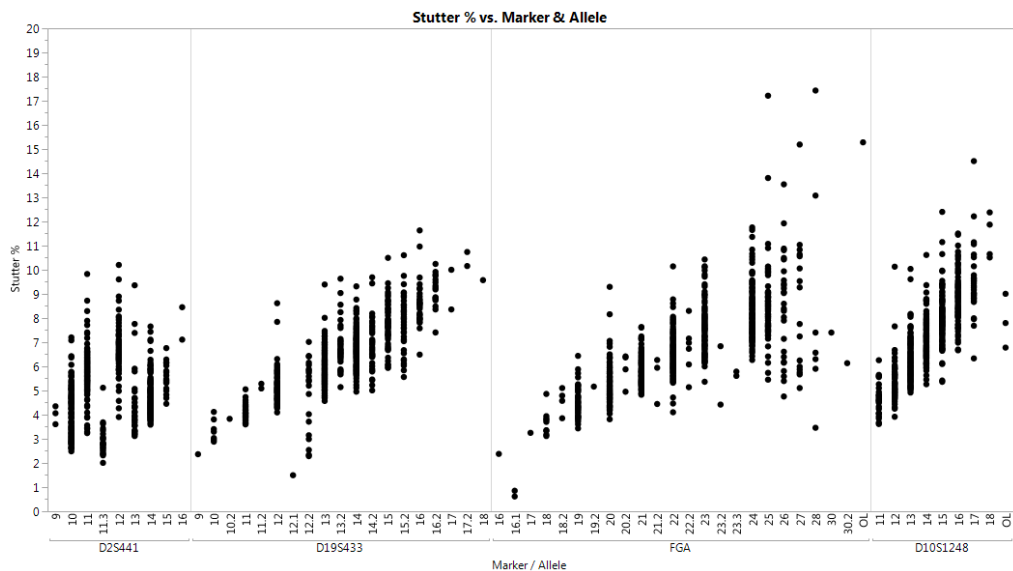


Figure 25 Stutter percentages for the Verifiler™ Plus kit TED™ dye (yellow) loci: D2S441, D19S433, FGA, and D10S1248.

Note: In the electropherogram plot, yellow is displayed as black.

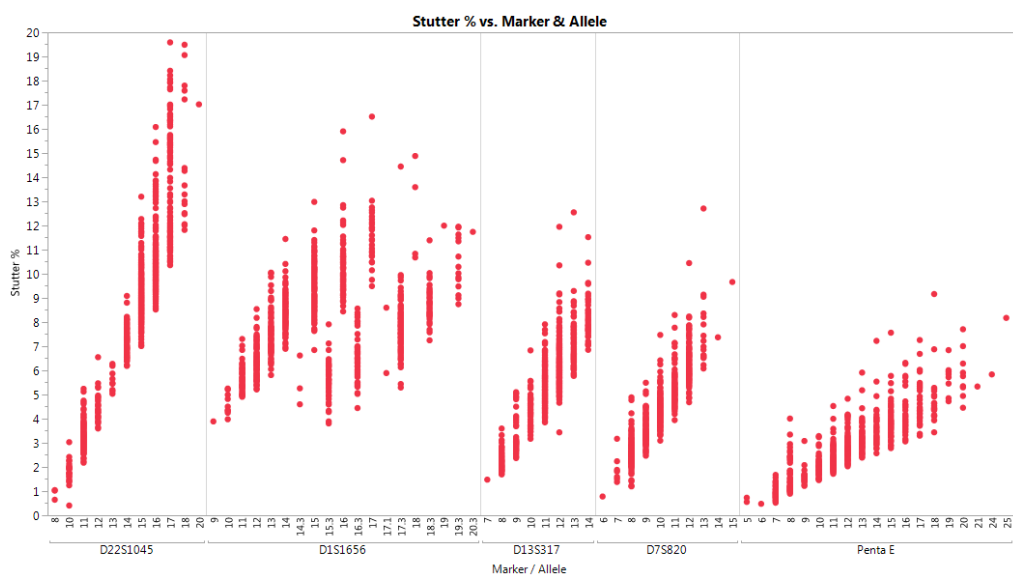


Figure 26 Stutter percentages for the Verifiler™ Plus kit TAZ™ dye (red) loci: D22S1045, D1S1656, D13S317, D7S820, and Penta E.

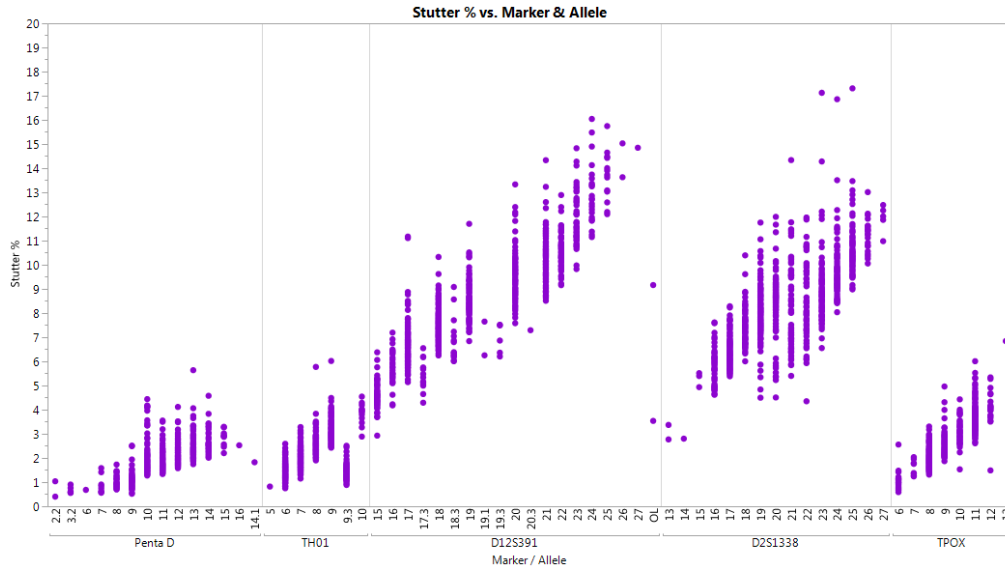


Figure 27 Stutter percentages for the Verifiler™ Plus kit SID™ dye (purple) loci: Penta D, TH01, D12S391, D2S1338, and TPOX.

## Plus-stutter peaks at the D22S1045 locus

The D22S1045 STR locus in the Verifiler™ Plus kit is a trinucleotide repeat locus, and shows an elevated level of plus stutter (Figure 28). Other loci, such as FGA, may also exhibit elevated plus stutter.

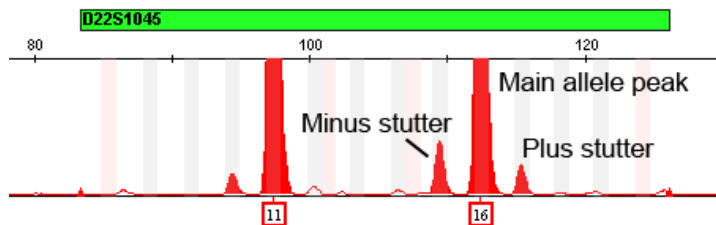


Figure 28 Verifiler™ Plus kit electropherogram showing minus and plus stutter associated with the D22S1045 STR locus. Data produced on a 3500xL Genetic Analyzer.

STR loci such as D1S1656 (Figure 29) 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™ ID-X Software for analysis of Verifiler™ Plus kit data contains a minus 2-bp stutter filter for D1S1656, as well as filters for commonly observed plus stutter, to prevent these peaks from being called in normal profiles.

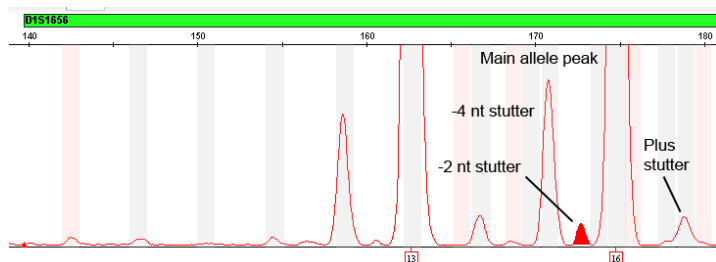


Figure 29 Example of non-standard stutter peaks in the D1S1656 locus. In addition to the main allele peaks and standard (–4-bp) stutter peaks, –2-bp minus stutter peaks and +4-bp plus stutter peaks can also be observed. Data produced on a 3500xL Genetic Analyzer.



## Stutter filter settings provided with the GeneMapper™ ID-X Software

The stutter filter settings are shown in Table 8. The data used to derive the settings are shown in Figure 23 through Figure 27. 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 minimized the occurrence of non-filtered stutter peaks while allowing maximum peak-detection sensitivity was determined.

---

**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 8 Percentages used in the stutter filters included with the GeneMapper™ ID-X Software**

**Note:** Penta D and Penta E do not require plus stutter filters because their stutter levels are so low.

Locus	Minus stutter (%)	Plus stutter (%)
D3S1358	12.61	3.74
vWA	11.86	4.46
D16S539	11.19	4.05
CSF1PO	11.00	5.42
D6S1043	12.05	4.58
D8S1179	11.54	4.85
D21S11	13.83	5.04
D18S51	13.73	5.96
D5S818	10.90	4.64
D2S441	9.71	3.71
D19S433	11.00	5.91
FGA	14.01	7.96
D10S1248	12.48	3.13
D22S1045	19.04	8.63
D1S1656	15.20	6.40
D13S317	11.00	6.25
D7S820	9.83	4.27
Penta E	7.99	—

**Table 8 Percentages used in the stutter filters included with the GeneMapper ID-X Software (continued)**

Locus	Minus stutter (%)	Plus stutter (%)
Penta D	4.59	—
TH01	5.55	3.95
D12S391	15.61	6.10
D2S1338	16.69	8.00
TPOX	5.82	3.49

## 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 observation

The efficiency of +A addition is related to the particular sequence of the DNA at the 3' end of the PCR product. The 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 for 5 minutes at 60°C.

This final extension step gives the DNA polymerase additional time to complete +A addition to all double-stranded PCR products. Examples of incomplete and normal +A addition are shown in Figure 30. 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.

If the amount of input DNA is greater than the 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. Other adverse conditions may also result in incomplete adenylation, for example: PCR inhibition, expired PCR reagents, or not completing the final hold step during thermal cycling.

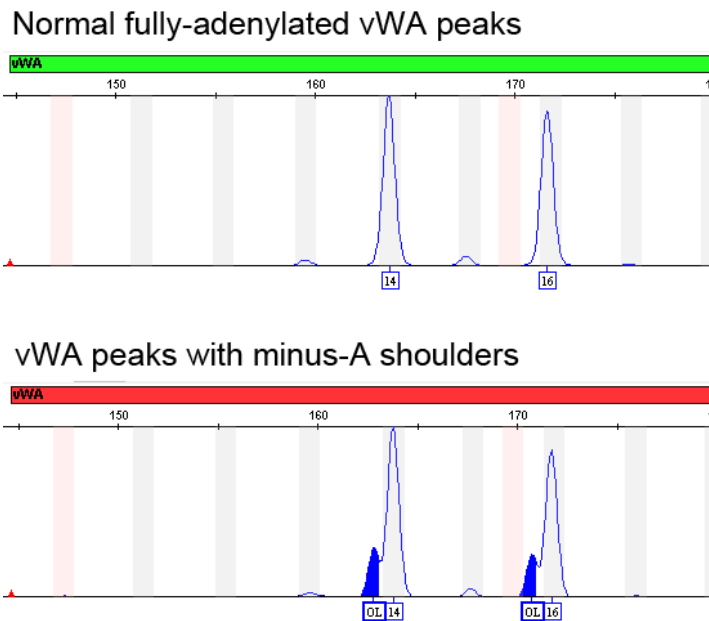


Figure 30 The top panel shows normal vWA peaks that are fully adenylation, with no visible shouldering and correct genotype calls. The bottom panel shows vWA peaks in which adenylation was incomplete, resulting in minus-A shoulders that were called as off-ladder (OL) peaks by the GeneMapper™ *ID-X* Software.

## Extra peaks: Artifacts

### Artifact definition

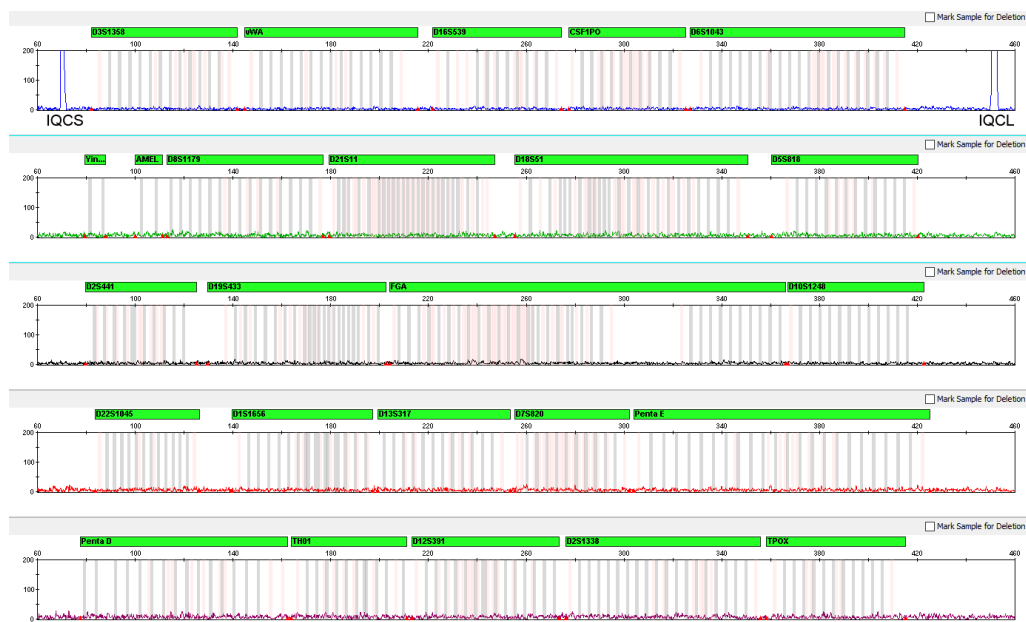
Artifacts and anomalies are seen in all molecular biological systems. Artifacts are typically reproducible. Anomalies are non-reproducible, intermittent occurrences that are not observed consistently in a system (for example, spikes and baseline noise).

### Dye artifact observation

Because of improvements in PCR primer manufacturing processes, the incidence of artifacts has been greatly reduced in the Verifiler™ Plus kit. Internal population studies show that kit electropherograms are free of reproducible dye artifacts in the kit read region of 64–458 bp. Two exceptions are as follows:

- A low-level 113–117-bp dye artifact in the VIC™ dye channel that has been detected.
- A low-level ~66-bp 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.

The low baseline-level fluorescence that is observed in a typical negative control PCR is shown in Figure 31. However, it is important to consider noise and other amplification-related artifacts when interpreting data.



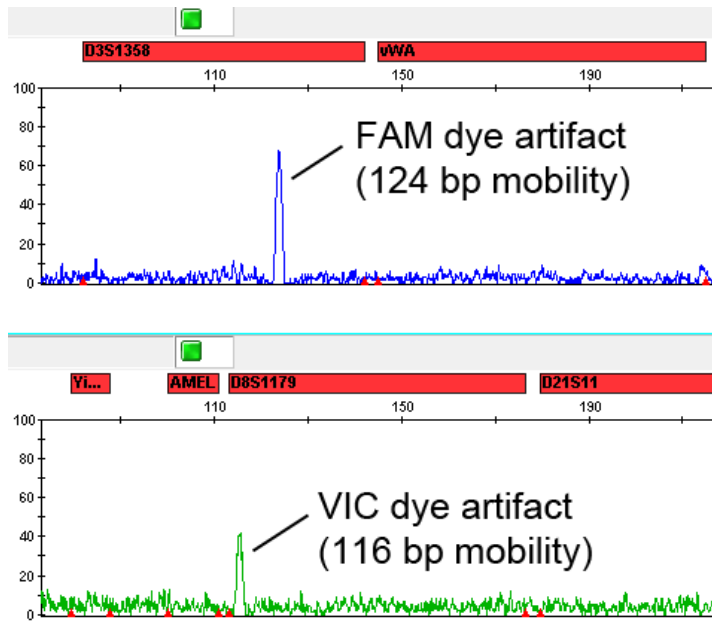
**Figure 31** Examples of fluorescence background in data produced on a 3500xL Genetic Analyzer (Y-axis scale 0–200 RFU) (updated formulation)

Note the two FAM™ dye-labeled IQC peaks (IQCS and IQCL) that should normally be present in negative samples.

Some small PCR artifacts were occasionally observed in negative control reactions prepared with the Verifiler™ Plus kit reagents that were subjected to long-term storage at elevated temperature as part of stability testing. The most prominent and consistent artifacts that were sometimes observed were:

- An artifact sizing at 120–130 bp in the FAM™ dye channel
- An artifact sizing at 110–120 bp in the VIC™ dye channel

For examples of these dye artifacts, see Figure 32. Such artifacts may form spontaneously upon long-term storage of PCR reagents, but their formation can be minimized by following the recommended storage conditions.



**Figure 32** Examples of FAM™ dye and VIC™ dye artifacts in negative (NTC) reactions with Verifiler™ Plus kit reagents that were subjected to long-term storage at elevated temperature  
In this example, both artifact peaks are below 100 RFU.

## Characterization of loci

### SWGDM guideline 3.1

“The basic characteristics of a genetic marker should be determined and documented.” (SWGDM, December 2016)

### Loci in this kit

This section describes basic characteristics of the 23 autosomal STR loci, Y indel locus, and sex-determining marker (Amelogenin), that are amplified with the Verifiler™ Plus 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 102-bp and 108-bp 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 81-bp or a 87-bp 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.) Most of the STR loci present in the Verifiler™ Plus kit contain tetranucleotide repeat units. However, the kit also contains one locus with trinucleotide repeats (D22S1045) and two loci with pentanucleotide repeats (Penta D and Penta E). 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 Verifiler™ Plus 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 Verifiler™ Plus 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

Four sets of STR loci in the Verifiler™ Plus kit are located on the same chromosome arms, as shown in Table 9. Of these, the most closely spaced are vWA and D12S391, which are located 6.3 million bp apart on the p arm of chromosome 12. Linkage disequilibrium analysis was conducted on vWA and D12S391 genotype results from 1,034 individuals of three ethnic groups (350 African-American, 349 Caucasian, and 335 Hispanic) using the Linkage Disequilibrium module of GenePop software version 4.0.10 (Raymond and Rousset, 1995; Rousset, 2008).

The results of this analysis (not shown) indicated that there is no statistically significant linkage disequilibrium found between vWA and D12S391.

However, an inheritance 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 (Budowle, *et al.*, 2010).

For examples of additional research on this topic, see: Burri, *et al.*, 2015; Liu, *et al.*, 2013.

**Table 9** STR loci in the Verifiler™ Plus kit located on the same chromosome arms

Locus	Chromosome location <sup>[1]</sup>		Distance apart (Mb)
	Map	Mb units	
D5S818	5q23.2	123.139	26.297
CSF	5q33.1	149.436	
vWA	12p13.31	5.963	6.378

Table 9 STR loci in the Verifiler Plus kit located on the same chromosome arms (continued)

Locus	Chromosome location <sup>[1]</sup>		Distance apart (Mb)
	Map	Mb units	
D12S391	12p13.2	12.341	6.378
D21S11	21q21.1	19.476	24.404
Penta D	21q22.3	43.88	
TPOX	2p25.3	1.472	66.741
D2S441	2p14	68.213	

<sup>[1]</sup> STR locus mapping data was obtained from the NCBI Map Viewer [http://www.ncbi.nlm.nih.gov/projects/mapview/map\\_search.cgi?taxid=9606](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) >0.05 were considered to indicate that linkage disequilibrium between the loci within the population tested was not statistically significant.

## Species specificity

### SWGAM 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.” (SWGAM, December 2016)

### Nonhuman study observation

The Verifiler™ Plus kit provides the required specificity for detecting human alleles. Species specificity testing was performed to show 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 kit:

- **Primates**—Chimpanzee, orangutan, and macaque (0.5 ng each)
- **Non-primates**—Mouse, dog, sheep, pig, rabbit, cat, horse, rat, and cow (5.0 ng each)
- **Microorganisms**—*Streptococcus salivarius*, *Neisseria gonorrhoea*, *Bacillus subtilis*, *Candida albicans*, *Lactobacillus delbrueckii*, *Escherichia coli*, and *Staphylococcus aureus* (pooled genomic DNAs, with approximately 100,000 copies of DNA from each species, per reaction)

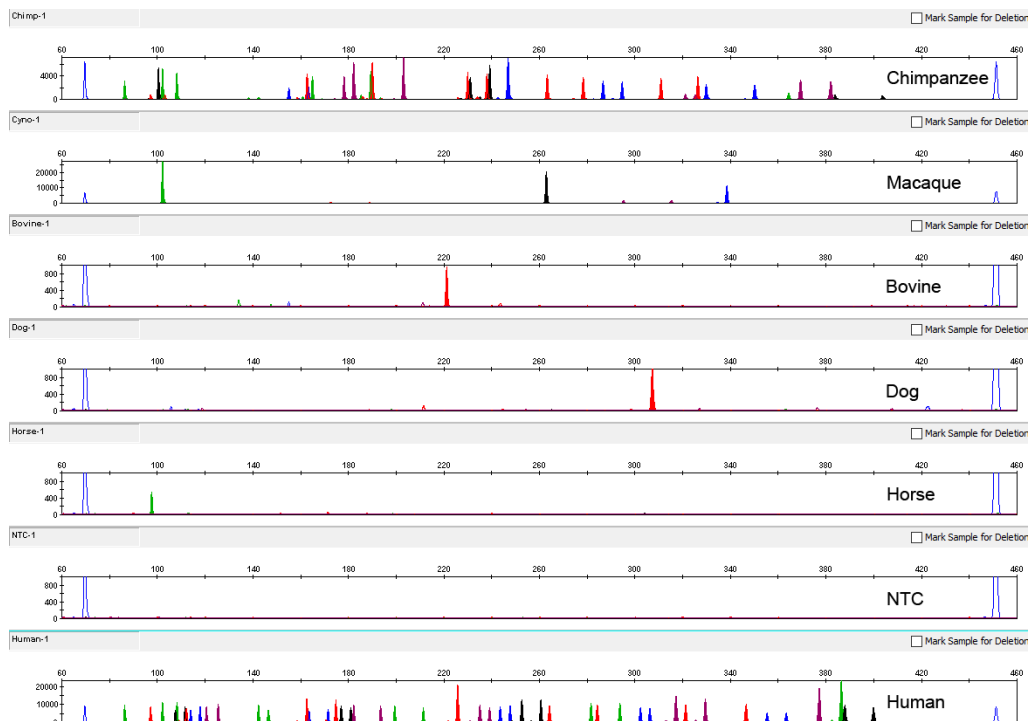
Results were evaluated for the presence of any amplified peaks that would indicate cross-reactivity of the kit with any of these non-human species.

Primate species (chimpanzee, macaque) produced more extensive cross-reactivity than non-primate species, and for a closely related species like chimpanzee, many peaks fell into allele bins.

A few of the more distantly related mammals (cow, dog, horse, hamster, and sheep) produced peaks over the 175 RFU threshold. No peaks fell into human STR locus bins, and would therefore not be confused with human STR alleles. None of the remaining species that were tested gave peaks over the detection threshold. These data show that the likelihood of obtaining an allelic profile from non-primates or microorganisms that is consistent with an allelic profile from a human sample is extremely low.

Example electropherogram results from the species specificity tests are shown in Figure 33.

The most significant cross-reactive peaks that were observed among non-human, non-primate, genomic DNAs (that is, peaks <175 RFU Peak Amplitude Threshold on the 3500xL Genetic Analyzer), with the original and updated formulations of the Verifiler™ Plus kit are provided in Table 10.



**Figure 33** Representative electropherograms for some species tested in a species specificity study. Data produced on a 3500xL Genetic Analyzer (updated kit formulation).

**Note:** The individual panes are at different magnifications (zoom function) on the Y-axis.



**Table 10** Observed cross-reactive peaks for non-human, non-primate animals for the original and updated formulations of the Verifiler™ Plus kit

Species	Dye channel	Size	Verifiler™ Plus kit—Original formulation		Verifiler™ Plus kit—Updated formulation	
			Average peak height	Genotype call	Average peak height	Genotype call
Dog	Blue	341 bp	214 RFU	D6 OL	—	—
	Red	307 bp	692 RFU	Penta E OL	716 RFU	Penta E OL
Horse	Green	98 bp	630 RFU	Out of Marker Range (OMR)	828 RFU	OMR
Hamster	Green	130 bp	243 RFU	D8 OL	—	—
Cow	Red	221 bp	399 RFU	D13 OL	489 RFU	D13 OL
Sheep	Purple	211 bp	345 RFU	TH01 OL	301 RFU	TH01 OL

## Sensitivity

### SWGAM 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.” (SWGAM, December 2016)

### Sensitivity observation

The recommended amount of input DNA for the Verifiler™ Plus kit is 0.5 ng for 29-cycle PCR, based on real-time PCR quantification with kits such as the Quantifiler™ Trio DNA Quantification Kit or the Quantifiler™ HP DNA Quantification Kit.

To determine the optimum input DNA amount, 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.

The amplification results of serially diluted DNA Control 007 (1.0–0.016 ng) are shown in Figure 34. Full profiles (46 human alleles) were consistently obtained at 0.125 ng, but occasional partial profiles resulted at lower concentrations.



Figure 34 Electropherograms for amplifications using 1,000, 500, 250, 125, 63, 31, and 16 pg of DNA Control 007 (updated formulation). Electrophoresis was performed on a 3500xL Genetic Analyzer. Note that because the DNA input is serially diluted by 2-fold, the Y-axis scale is also magnified by 2-fold for the smaller input amounts of DNA.

## Stability

### SWGAM 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.” (SWGAM, December 2016)

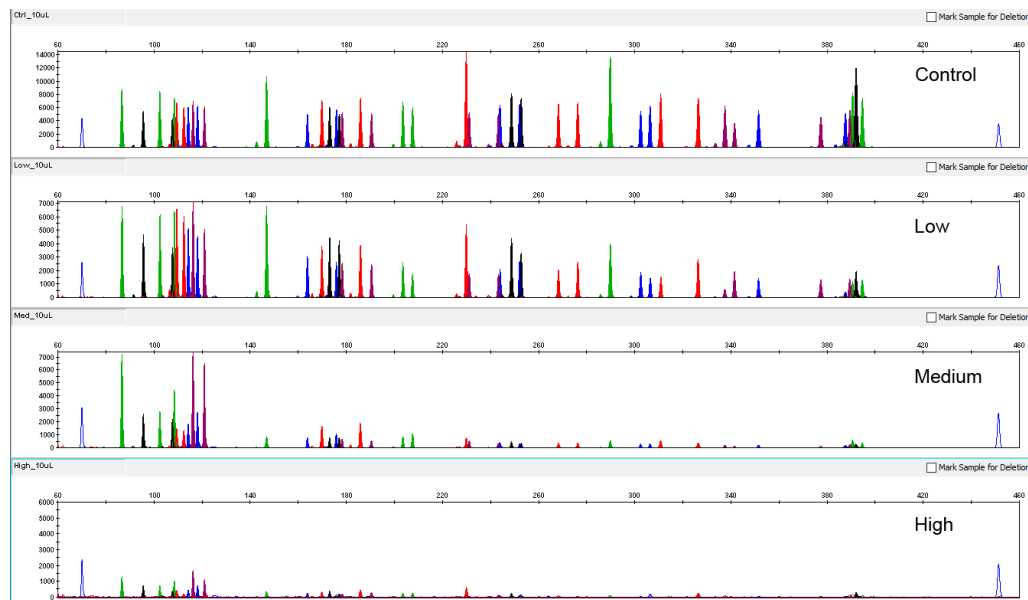
### Lack of amplification of some loci

As with any multi-locus system, the possibility exists that not every locus amplifies. This possibility is most often observed when the DNA sample contains PCR inhibitors or when the DNA sample has been severely degraded. Valuable information can be obtained from partial profiles.

## Degraded DNA

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

Degraded DNA was prepared to examine the potential for preferential amplification of loci. High-molecular-weight DNA was sonicated and incubated with increasing doses of DNase I (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 Verifiler™ Plus 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.



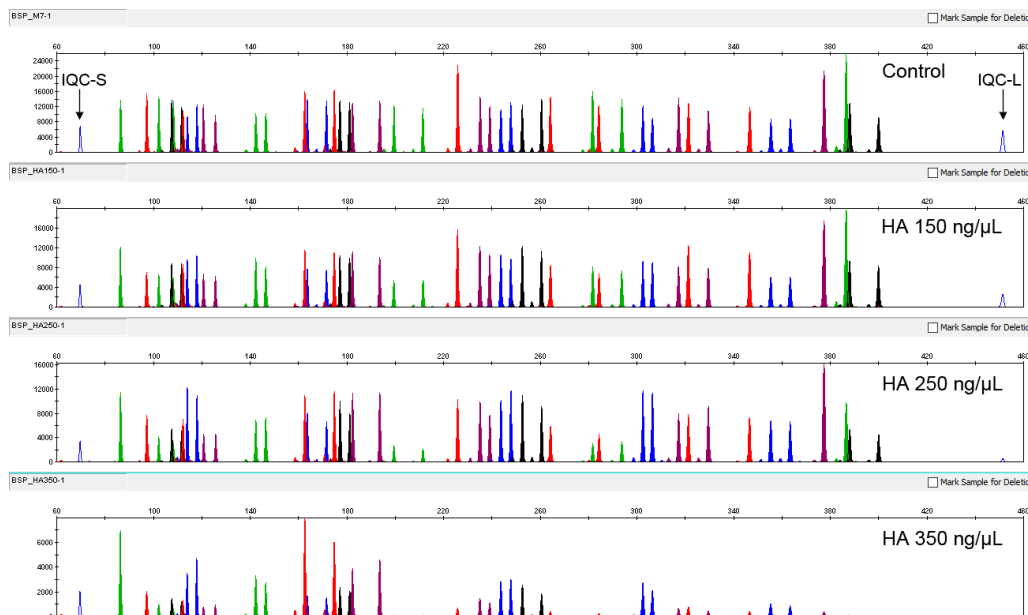
**Figure 35 Amplification of a single donor DNA sample sonicated and incubated with increasing doses of DNase I (updated formulation)**

The panels show the control (undegraded), low-, medium-, and high-degraded samples. Note that the Y-axis scale is magnified for more degraded samples, which generate lower peak heights.

## Effect of inhibitors

Because of the substrates and environments from which they are collected, forensic DNA samples may be contaminated with substances that inhibit PCR amplification. The PCR chemistry for the Verifiler™ Plus kit was optimized to be as robust as possible in the presence of such inhibitors. Model inhibitors tested during developmental validation were: hematin, a representative compound often extracted from blood stains (DeFranchis *et al.*, 1988; Alkane *et al.*, 1994); humic acid, a simulant for soil-based inhibition; and tannic acid, a compound often co-extracted from leather sample substrates. 0.5 ng of DNA Control 007 was amplified in the presence of increasing concentrations of each model inhibitor. All samples were then amplified using the standard 29-cycle "touchdown" thermal cycling program.

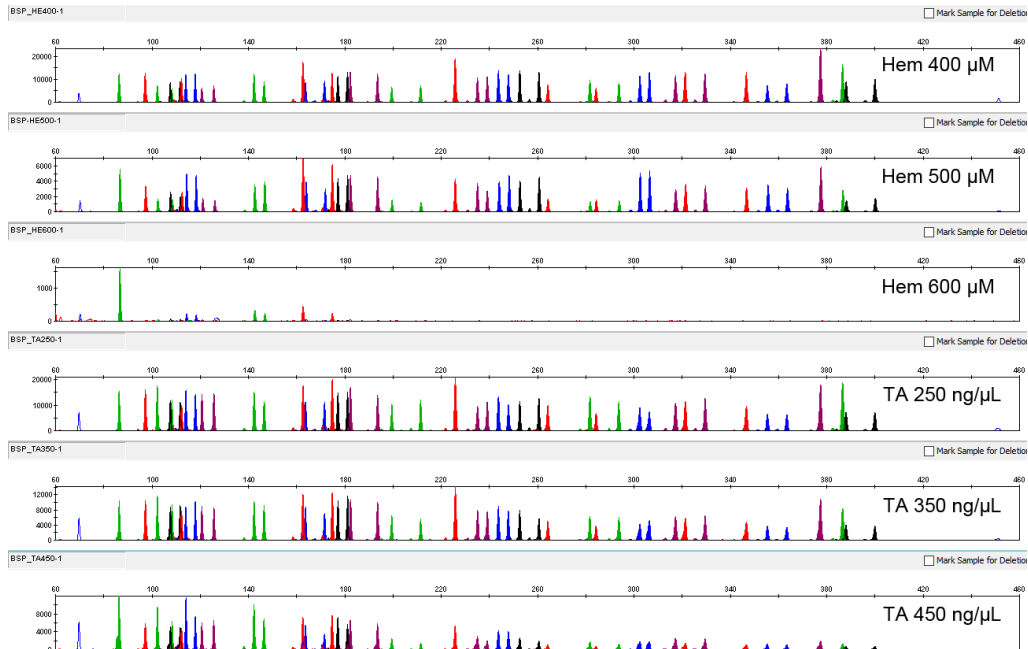
Representative electropherograms from the PCR inhibitor studies with humic acid, hematin, and tannic acid are shown in Figure 36 and Figure 37. Inhibitor final concentrations in 25- $\mu$ L PCR reactions and average allele counts obtained in replicate reactions with the Verifiler™ Plus kit are shown in Table 11.



**Figure 36** Electropherograms show the results of testing the Verifiler™ Plus kit with the humic acid PCR inhibitor (updated formulation)

The top panel is the experimental control with 0.5 ng (total) of DNA Control 007 without humic acid. The remaining panels show 0.5 ng of DNA Control 007 with increasing levels of humic acid (150, 250, and 350 ng/μL). Note the behavior of the IQCS and IQCL peaks in response to increasing levels of inhibition, with the IQCL peak decreasing in peak height and eventually disappearing as the humic acid concentration increases.

**Note:** The individual panes are at different magnifications (zoom function) on the Y-axis.



**Figure 37** Example electropherograms show the results of testing hematin and tannic acid PCR inhibitors (updated formulation)

Each sample contained 0.5 ng (total) of human male DNA Control 007 with varying levels of hematin or tannic acid. Note that the IQCL peak was significantly reduced or absent while the IQCS peak still remained in the inhibitor-containing samples.

**Note:** The individual panes are at different magnifications (zoom function) on the Y-axis.

**Table 11** Verifiler™ Plus kit performance with PCR inhibitors (n=4) (updated formulation)

Sample	Inhibitor concentration	Mean allele count
Control (no inhibitor)	0	46
Hematin low	400 μM	46
Hematin medium	500 μM	46
Hematin high	600 μM	8.75
Humic acid low	150 ng/μL	46
Humic acid medium	250 ng/μL	46
Humic acid high	350 ng/μL	42.5
Tannic acid low	250 ng/μL	46
Tannic acid medium	350 ng/μL	46
Tannic acid high	450 ng/μL	46

## Mixture studies

### SWGDM guideline 3.8

“The ability to obtain reliable results from mixed source samples should be determined.” (SWGDM, December 2016)

### Mixture study overview

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 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.

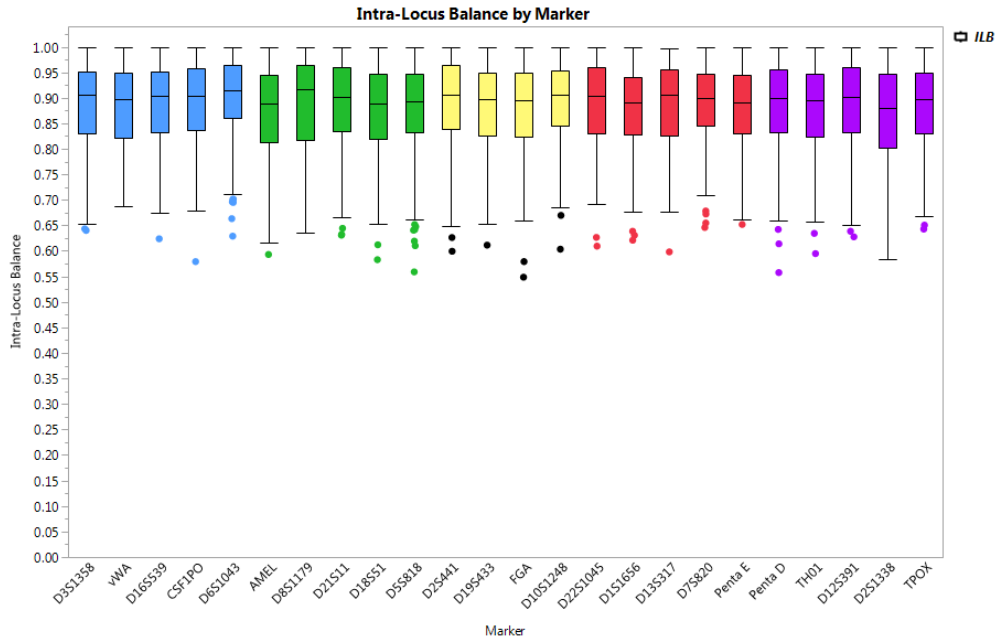
The peak height ratio is defined as the height of the lower peak (in RFU) divided by the height of the higher peak (in RFU), expressed as a percentage.

If an unusually low peak height ratio is observed for one locus, and there are no other indications that the sample is a mixture, re-amplify 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 Verifiler™ Plus kit loci in single-source human population database samples are shown in Figure 38. The population samples that are used are listed in “Population samples used in these studies” on page 99.



**Figure 38** Heterozygote ratios for ~0.5 ng of input DNA. Intra-locus peak height ratio is calculated per individual profile by dividing the lower peak by the higher peak in each heterozygote allele pair, per marker. Box plot boxes and data points are colored according to the dye channel. 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.

## Resolution of genotypes in mixed samples

A sample that contains DNA from two sources can comprise (at a single locus) any of the following seven genotype combinations:

- 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™ instruments provides additional valuable data to aid in resolving mixed genotypes.

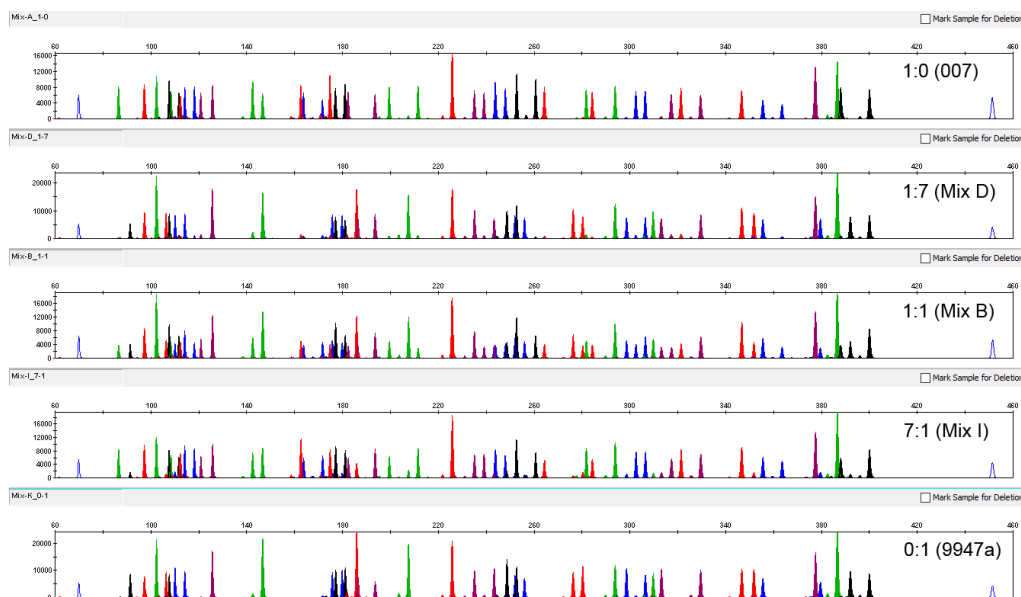
Ultimately, the likelihood that any sample is a mixture must be determined by the analyst in the context of each particular case, including the information provided from known reference samples.

## Limit of detection of the minor component

Mixtures of two DNA samples were examined at various M:F ratios (0:1, 1:1, 1:3, 1:7, 1:15, 1:30, 30:1, 15:1, 7:1, 3:1 and 1:0). The total amount of genomic input DNA mixed at each ratio was 0.5 ng. The samples were amplified in a ProFlex™ 96-well PCR System, then electrophoresed and detected using an 3500xL Genetic Analyzer.

The results of the mixed DNA samples are shown in Figure 39. The two human genomic DNAs, male DNA Control 007 and female Control DNA 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 3/3 replicates of both samples giving full profiles of the respective minor contributor. 15:1, 30:1, 1:30, and 1:15 ratios resulted in partial profiles for the minor component (Table 13). The genotypes of each contributor DNA are shown in Table 12.

Mixture sample compositions and mean STR allele counts from Verifiler™ Plus kit assays are shown in Table 13.



**Figure 39** Amplification of DNA mixtures at various ratios (updated formulation). Panels show electropherograms for 1:0 (male DNA Control 007 only), 1:7 mixture, 1:1 mixture, 7:1 mixture, and 0:1 (female Control DNA 9947A only).



**Table 12 Genotypes of mixed DNA samples (updated formulation). (Asterisks indicate alleles that are unique for each individual, and also apart from plus- or minus-stutter positions of other alleles present in the individuals.)**

Locus	Male 007 genotype		Female 9947A genotype	
	Allele 1	Allele 2	Allele 1	Allele 2
D3S1358	15	16*	14	15
vWA	14*	16	17*	18*
D16S539	9*	10	11*	12*
CSF1PO	11	12	10	12
TPOX	8	—	8	—
Y indel	2*	—	—	—
AMEL	X	Y*	X	—
D8S1179	12	13	13	—
D21S11	28*	31*	30	—
D18S51	12*	15	15	19*
D2S441	14	15*	10*	14
D19S433	14	15	14	15
TH01	7	9.3	8*	9.3
FGA	24	26*	23	24
D22S1045	11	16*	11	14*
D5S818	11	—	11	—
D13S317	11	—	11	—
D7S820	7*	12*	10*	11
D10S1248	12	15	13*	15
D1S1656	13*	16*	18.3*	—
D12S391	18	19	18	20*
D2S1338	20*	23	19	23
D6S1043	12	14*	12	18*
Penta E	7*	12	12	13*
Penta D	11	12	12	—

**Table 13 Mixture sample compositions and mean STR allele counts from Verifiler™ Plus kit assays (updated formulation)**

Mixture sample	Ratio (A:B)	DNA amounts (ng/reaction)		Mean allele count <sup>[1]</sup>	
		Male 007 ("A")	Female 9947A ("B")	Male 007 ("A")	Female 9947A ("B")
A	1:0	0.500	0.000	18.00	0.00
B	1:1	0.250	0.250	18.00	14.00
C	1:3	0.125	0.375	18.00	14.00
D	1:7	0.063	0.438	18.00	14.00
E	1:15	0.031	0.469	15.33	14.00
F	1:30	0.016	0.484	12.67	14.00
G	30:1	0.484	0.016	18.00	7.33
H	15:1	0.469	0.031	18.00	10.67
I	7:1	0.438	0.063	18.00	14.00
J	3:1	0.375	0.125	18.00	13.67
K	0:1	0.000	0.500	0.00	14.00

<sup>[1]</sup> Mixture sample compositions and STR allele counts from Verifiler™ Plus kit assays. The 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. The full profiles of unique alleles in the DNA Control 007 and Control DNA 9947A DNAs were 18 and 14, respectively.

## Population data

### SWGAM guideline 3.7

“The distribution of genetic markers in populations should be determined in relevant population groups.” (SWGAM, December 2016)

### Loci in the kit

The Verifiler™ Plus PCR Amplification Kit was designed and optimized primarily for the analysis of forensic casework samples. It was designed as a companion kit to the Verifiler™ Express PCR Amplification Kit (Cat. No. A32014) and the Huaxia™ Platinum™ PCR Amplification Kit (Cat. No. A31323), and contains the same set of 25 loci.

## Population distribution

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.

## Population samples used in these studies

The Verifiler™ Plus kit has high genotypic concordance (>99.5%) to its databasing counterparts, the Verifiler™ Express kit and Huaxia™ Platinum™ kit (all three kits contain the same 25 loci). An initial population study of individuals of different ethnic populations within the United States and China was conducted with 2,475 individuals. This initial study was the basis for the allele frequency tables and associated random match probability tables at the end of this section.

After the minor redesign of certain primers in the Verifiler™ Plus kit multiplex to reduce the incidence of artifact peaks in samples with high levels of bacterial DNA, a new population study was performed to check:

- Concordance between the original Verifiler™ Plus kit and the redesigned kit
- Concordance with the GlobalFiler™ kit reference genotypes

As with the original population study, the study with the updated formulation included in-house testing of domestic United States individuals as well as population samples collected and tested by laboratories in China. See Table 14.

**Table 14** Population samples (updated formulation)

Sample source (population)	Total samples	Sample type	No. of samples by ethnic origin
In-house testing of domestic United States individuals	1,308	<ul style="list-style-type: none"> <li>• Extracted DNA from 1,108 blood samples</li> <li>• 200 samples from blood cards that were tested in direct amplification mode</li> </ul>	<ul style="list-style-type: none"> <li>• African-American—285</li> <li>• Asian—316</li> <li>• Caucasian—322</li> <li>• Hispanic—385</li> </ul>
Chinese laboratories	987	Blood cards that were tested in direct amplification mode	Individuals of Chinese origin—987

## Concordance studies

Concordance between the original Verifiler™ Plus kit and the redesigned kit was 100% for 2,288 samples tested in-house and externally. When genotype results were compared between the Verifiler™ Plus kit and the GlobalFiler™ kit, the following level of concordance was observed:

Sample source (population)	Sample type	Level of discordance
In-house testing of domestic United States individuals	Extracted DNAs from 1,108 blood samples	1,099/1,108 (discordances observed in D10S1248 and D3S1358)
	200 samples from blood cards that were tested in direct amplification mode	200/200 (0 discordances observed)
Chinese laboratories	Blood cards that were tested in direct amplification mode	986/987 (1 discordance observed in D5S818)

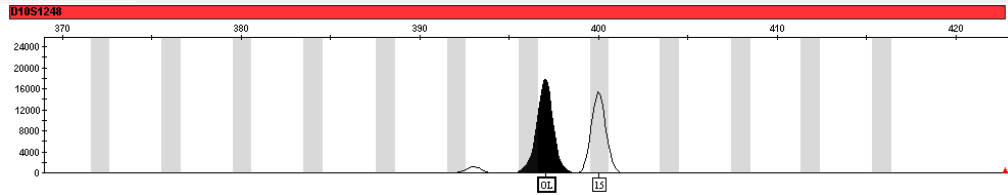
Among the in-house extracted samples, several instances of discordance were observed that were caused by a 1-bp insertion in the flanking region of D10S1248. Affected individuals genotyped as off-ladder (OL). The same individuals tested with the GlobalFiler™ kit did not give the OL genotypes. All individuals that contained the 1-bp insertion were of African descent, and subsequent database searches confirmed that this indel is found almost exclusively in individuals of African ancestry.

### D10S2148 in the Verifiler™ Plus kit

The Verifiler™ Express kit and Huaxia™ Platinum™ kit were designed for database workflows that involve pristine single-source samples. The Verifiler™ Plus kit was designed for casework workflows that involve challenging samples. Although the three kits are complementary, the Verifiler™ Plus kit was modified to optimize casework performance and discrimination power with degraded samples. In the Verifiler™ Plus kit, more discriminating markers (such as the Penta D marker) were given priority for amplification and located in the smaller molecular weight regions, while the D10S1248 marker, because of its lower discrimination power, was positioned in the larger molecular weight region.

To move the D10S1248 marker from a smaller base pair size in the SID™ dye channel to the larger base pair size in the TED™ dye channel, the flanking region that borders the primer binding and STR sites was extended. By sequencing additional base pairs in the flanking region of the D10S1248 marker, the opportunity to observe rare, sample-specific mutations increases.

In the case of six African-American (out of 330) and one Hispanic (out of 393) population samples, an insertion in the flanking region for the D10S1248 marker caused a 1-bp difference when comparing results from the Verifiler™ Plus kit to the Verifiler™ Express kit, Huaxia™ Platinum™ kit, and GlobalFiler™ kit (Figure 40). This was seen with alleles 13, 14, and 15 reported as 13.1, 14.1, and 15.1. The discrepancy has not been observed in any other population samples tested (1,335 total combined Caucasian and Asian samples). Subsequent sequencing of the sample showed that it is consistent with African lineage. This D10S1248 insertion is observed in the original and updated formulations of the Verifiler™ Plus kit.



**Figure 40** D10S1248 marker from an African American population sample amplified with the Verifiler™ Plus kit. The off-ladder (OL) peak sizes at the 14.1 base pair position because of a single base pair insertion in the flanking region between the primer and STR region. When amplified with the Verifiler™ Express kit, Huaxia™ Platinum™ kit, and GlobalFiler™ kit, the genotype at this marker would be 14, 15.

## Supporting data

Additional studies were performed at the National Institute of Standards and Technology (NIST, C.R. Steffan). Those studies confirmed the above conclusions, showing the 1-bp difference in nine out of 355 African-American samples.

For expanded frequency analysis of this insertion, we examined the sequencing data for the specific insertion at the D10S1248 marker from the following organizations:

- Broad Institute ([https://gnomad.broadinstitute.org/variant/10-131092326-T-TA?dataset=gnomad\\_r2\\_1](https://gnomad.broadinstitute.org/variant/10-131092326-T-TA?dataset=gnomad_r2_1))
- 1,000 Genomes Project (<https://www.internationalgenome.org/about>)
- NCBI ([https://www.ncbi.nlm.nih.gov/snp/rs567190580#frequency\\_tab](https://www.ncbi.nlm.nih.gov/snp/rs567190580#frequency_tab))

Out of 5,044 total African-related samples, there were 80 observations. Out of 771 total Hispanic-related samples, the insertion was not observed.

## Next steps and considerations

The potential to observe discrepant genotypes among STR PCR kits that use different primer sequences is a known phenomenon in the forensic community. Depending on workflow, interpretation protocols, and database search parameters, each laboratory may take a different approach to handling discrepancies across STR PCR kits.

If your laboratory has an established protocol for handling these discrepancies, we recommend that you follow that protocol for this occurrence. If your laboratory does not have an established protocol, we recommend that you consider different analysis and database search approaches, some of which include:

- Evaluate previously generated Verifiler™ Plus kit files that contain 0.1 microvariant calls at the D10S1248 marker.
- Confirm the genotype through amplification with a secondary STR PCR kit, such as the GlobalFiler™ kit.
- Exclude the D10S1248 marker from upload to a database if the 0.1 microvariant is observed.
- Perform database searches that include one mismatch when a database is composed of profiles from different STR PCR kits.

## Probability of identity

### Probability of identity definition

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

### Probability of identity observation

The autosomal STR allele frequencies at Verifiler™ Plus kit loci by population group are shown in Figure 41 through Figure 45.

---

**Note:** These tables show observed allele frequencies among different population groups. To ensure a conservative approach when performing statistical calculations such as profile frequency estimates, minimum allele frequencies should be used for very rare alleles that may be insufficiently sampled in a population. For autosomal markers, minimum allele frequencies are typically calculated as  $5/2N$ , where  $N$  is the number of individuals sampled in a given population (Butler, 2010).

---

The Random Match Probability (RMP) values of the Verifiler™ Plus kit loci (individually and combined) are shown in Table 15.

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA	
Sample Size	2463	2462	2455	2463	2461	2462	2460	2463	2462	2462	2447	2463	2458	2458	2463	2463	2463	2461	2462	2460	2454	2438	2463	
4							0.0002																	
5							0.0002												0.0002	0.0467	0.0002			
6	0.0002						0.0002									0.0002			0.0039	0.0004	0.1057	0.0002		
7	0.0037			0.0008	0.0002		0.0002							0.0173	0.0008	0.0032			0.0104	0.0124	0.2675	0.0016		
8	0.0018	0.0022		0.2542	0.0128			0.0006						0.0028	0.0035	0.1592	0.0022		0.0471	0.0053	0.0607	0.5455		
9	0.0477			0.1326	0.2485		0.0002					0.0006		0.0663	0.0012	0.0666	0.0022		0.3148	0.0116	0.4817	0.1081		
9.1												0.0199				0.0035								
9.2																0.0008								
9.3																					0.0591			
10	0.2326	0.0010		0.1435	0.1117	0.0012	0.0002	0.0006		0.0002		0.2485		0.1937	0.0298	0.1661	0.1139		0.1265	0.0488	0.0247	0.0277	0.0004	
10.1												0.0014												
10.2							0.0004																	
10.3																							0.0002	
11	0.2420	0.0028		0.2426	0.2802	0.0043	0.0028	0.0607		0.2279		0.3388	0.0002	0.3216	0.1407	0.3244	0.0751		0.1877	0.1163	0.0002	0.2859	0.0002	
11.1																0.0002								
11.2							0.0004																	
11.3												0.0406												
12	0.3871	0.0701		0.1657	0.2263	0.0426	0.0431	0.0585		0.0032		0.1665	0.0012	0.2496	0.1360	0.2328	0.1242		0.1446	0.1195	0.0002	0.0287	0.0002	
12.1	0.0002						0.0002																	
12.2							0.0041																	
12.3												0.0012			0.0006									
13	0.0696	0.3400		0.0491	0.1052	0.2074	0.2677	0.1076		0.0051	0.0004	0.0264	0.0024	0.1391	0.1240	0.0384	0.2288	0.0002	0.1188	0.0543		0.0014	0.0008	
13.2							0.0402																	
13.3												0.0004												
14	0.0132	0.2400	0.0002	0.0110	0.0144	0.1881	0.2500	0.0818		0.0268		0.1447	0.0391	0.0077	0.1236	0.0041	0.1908	0.0004	0.0329	0.0876		0.0004	0.2245	
14.2							0.1037																	
14.3														0.0002										
15	0.0018	0.2226	0.0155	0.0004	0.0006	0.1673	0.0858	0.2722		0.2685		0.0104	0.3519	0.0014	0.0148	0.0002	0.1724		0.0097	0.1035			0.0335	

Figure 41 Combined China populations (minimum allele frequency = 0.0010)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA	
15.2							0.1370																0.0002	
15.3								0.0047																
16		0.0981	0.0094	0.0002		0.1355	0.0189	0.2259		0.2547	0.0119	0.0006	0.3255		0.0032		0.0771	0.0006	0.0032	0.0785		0.0002	0.1908	
16.1								0.0002																
16.2								0.0380												0.0002				
16.3								0.0085																
17		0.0213	0.1061			0.0810	0.0018	0.0855		0.1879	0.0719		0.2040		0.0376		0.0118	0.0008	0.0002	0.0772				0.2540
17.2							0.0039											0.0002						
17.3			0.0010					0.0583							0.0010									
17.4																				0.0002				
18		0.0018	0.2399			0.0437	0.0002	0.0108		0.0242	0.1030		0.0710		0.1726		0.0012	0.0254		0.0791				0.1841
18.1															0.0002									
18.2			0.0002				0.0006								0.0008									
18.3			0.0006					0.0191																
18.4																				0.0012				
19			0.1980			0.0469		0.0010		0.0016	0.1841		0.0045	0.0002	0.1439		0.0002	0.0453		0.0547				0.0962
19.2																		0.0002						
19.3			0.0004					0.0037							0.0002									
19.4																				0.0012				
20			0.1664			0.0303		0.0002			0.1165		0.0002		0.0540			0.0498		0.0437				0.0132
20.2																		0.0006						
20.3			0.0002					0.0002							0.0006									
21			0.1018			0.0217					0.0345				0.0071			0.1121		0.0209				0.0018
21.1						0.0002																		
21.2																		0.0035						
21.3															0.0024									
22			0.0843			0.0158					0.0499				0.0010			0.1772		0.0211				
22.2																		0.0049						
22.3			0.0002												0.0002									
23			0.0446			0.0077					0.2058							0.2117		0.0089				

Combined China populations (continued)



Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
23.2																		0.0102					
23.3											0.0002												
24			0.0187			0.0041			0.0002		0.1385							0.1784		0.0047			
24.2																		0.0069					
25			0.0094			0.0014					0.0666							0.1065		0.0016			
25.2																		0.0047					
26			0.0029			0.0006					0.0110							0.0427		0.0004			
26.2																		0.0010					
27			0.0002			0.0002			0.0028		0.0043							0.0116					
27.2																		0.0006					
28									0.0520		0.0012							0.0037					
28.2									0.0120														
29									0.2626		0.0002							0.0010					
29.2									0.0020														
30									0.2571														
30.1									0.0002														
30.2									0.0185														
30.3									0.0035														
31									0.0952														
31.2									0.0796														
31.3									0.0002														
32									0.0268														
32.2									0.1320														
33									0.0037														
33.2									0.0455														
34									0.0016														
34.2									0.0037														
35.2									0.0006														
39									0.0002														

Combined China populations (*continued*)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
Sample Size	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304	304
2.2																			0.1151				
3.2																			0.0263				
5																			0.0576	0.0789	0.0049		
6																0.0033			0.0115		0.1628	0.0822	
6.1																					0.0016		
7	0.0345													0.0033		0.0049			0.0230	0.1168	0.3618	0.0164	
7.3	0.0016																						
8	0.0905			0.0263	0.0329					0.0049		0.0016		0.0691		0.2039	0.0033		0.1003	0.1842	0.2023	0.3618	
9	0.0378	0.0016		0.0247	0.2155		0.0016	0.0016					0.0033	0.0214		0.1234	0.0033		0.1497	0.0493	0.1645	0.1891	
9.3																					0.0872		
10	0.2582			0.0280	0.1036	0.0033	0.0115	0.0132		0.0444		0.0855		0.0707	0.0148	0.3059	0.0362		0.1201	0.0395	0.0148	0.1020	
10.2							0.0016																
11	0.2253	0.0378		0.2878	0.2993	0.0049	0.0970	0.0510		0.1546		0.3536		0.2451	0.1053	0.2007	0.0609		0.1595	0.0658		0.2138	0.0033
11.2							0.0033																
11.3												0.0345											
12	0.2928	0.1414		0.4391	0.2007	0.0691	0.1086	0.0855		0.0658		0.1974		0.3438	0.2155	0.1332	0.1151		0.1266	0.1217		0.0329	
12.2							0.0395												0.0016				
13	0.0477	0.2204		0.1431	0.1332	0.0378	0.2681	0.1151		0.0016	0.0016	0.0378	0.0066	0.2270	0.0839	0.0230	0.1891		0.0789	0.1036		0.0016	0.0099
13.2						0.0033	0.0559																
14	0.0115	0.2829		0.0493	0.0148	0.0526	0.1941	0.2484		0.0806		0.2697	0.0872	0.0164	0.0789	0.0016	0.3421		0.0247	0.0691			0.0740
14.2						0.0033	0.0559																
14.3								0.0099															
15		0.1941	0.0691	0.0016		0.1628	0.0674	0.1776		0.2270	0.0033	0.0197	0.2780	0.0033	0.0444		0.1826		0.0049	0.0757			0.2253
15.1			0.0016																				
15.2							0.0444					0.0033											
15.3								0.0164															
16		0.0954	0.0526			0.1842	0.0148	0.0938		0.2039	0.0526		0.3322		0.0378		0.0559			0.0477			0.2615
16.1			0.0016															0.0033					

Figure 42 United States—African-American population (minimum allele frequency = 0.0082)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
16.2							0.0329																
16.3								0.0691															
17		0.0230	0.1678			0.1645		0.0280		0.1957	0.0921		0.2237		0.0905		0.0115			0.0280			0.1924
17.1			0.0016																				
17.2							0.0033																
17.3			0.0049					0.0609															
18		0.0033	0.2632			0.1447		0.0049		0.0214	0.0477		0.0609		0.1365			0.0099		0.0164			0.1398
18.2																		0.0082					
18.3			0.0082					0.0164															
19			0.1365			0.0938		0.0016			0.1563		0.0049		0.1118			0.0658		0.0016			0.0658
19.1			0.0049																				
19.2																		0.0033					
19.3			0.0033					0.0049															
20			0.1184			0.0461					0.1036				0.0625			0.0674		0.0016			0.0214
20.2						0.0016												0.0033					
20.3								0.0016															
21			0.0658			0.0197					0.1168				0.0049			0.1234					0.0066
21.2																		0.0016					
21.3			0.0016												0.0016								
22			0.0461			0.0049					0.1283				0.0016			0.1760					
22.2																		0.0016					
23			0.0362			0.0033					0.0987				0.0066			0.1612					
23.3																		0.0033					
24			0.0066								0.0938				0.0033			0.1908					
25			0.0099								0.0707							0.0938					
26											0.0280							0.0428					
27									0.0576		0.0066							0.0296					
28									0.2549									0.0115					
29									0.1579														
29.3									0.0016														

United States—African-American population (continued)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
30									0.2138									0.0016					
30.2									0.0132									0.0016					
31									0.0872														
31.2									0.0428														
32									0.0115														
32.2									0.0724														
33									0.0082														
33.2									0.0345														
34									0.0016														
35									0.0345														
36									0.0082														

United States—African-American population (*continued*)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
Sample Size	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376	376
5																				0.0426			
6																			0.0053	0.0013	0.1117		
7	0.0053			0.0013										0.0146	0.0040	0.0027			0.0080	0.0133	0.2739		
8	0.0013	0.0013		0.2301	0.0066			0.0013						0.0013	0.0027	0.1556	0.0027		0.0372	0.0027	0.0505	0.5133	
9	0.0545			0.1423	0.2779							0.0013		0.0572	0.0013	0.0505	0.0013		0.2979	0.0053	0.4973	0.1197	
9.1												0.0253				0.0027							
9.3																					0.0465		
10	0.2181	0.0013		0.1303	0.1184							0.2646		0.1915	0.0319	0.1569	0.0851		0.1423	0.0638	0.0186	0.0146	
10.1												0.0013											
11	0.2367			0.2766	0.2926	0.0040	0.0040	0.0598		0.2380		0.3511		0.3364	0.1316	0.3590	0.0718		0.1742	0.1104	0.0013	0.3245	
11.3												0.0372											
12	0.4122	0.0691		0.1689	0.1862	0.0346	0.0439	0.0465		0.0066		0.1423	0.0027	0.2380	0.1476	0.2327	0.1210		0.1569	0.1144		0.0253	
12.1	0.0013																						
12.2							0.0053																
12.3												0.0027											
13	0.0545	0.3471		0.0426	0.0997	0.2035	0.2620	0.0997		0.0040		0.0239		0.1516	0.1184	0.0372	0.2194	0.0013	0.1330	0.0585			
13.2							0.0572																
14	0.0093	0.2633		0.0080	0.0186	0.2074	0.2726	0.0811		0.0279		0.1370	0.0412	0.0080	0.1303	0.0027	0.1915		0.0306	0.1051		0.0027	0.2447
14.2							0.1170																
14.3														0.0013									
15	0.0066	0.2114	0.0160			0.1715	0.0864	0.3112		0.2660		0.0133	0.3684		0.0160		0.2008		0.0093	0.1077			0.0412
15.2							0.1024																
15.3								0.0040															
16		0.0864	0.0106			0.1316	0.0120	0.2101		0.2340	0.0080		0.3085		0.0066		0.0931		0.0053	0.0838			0.1662
16.2							0.0319																
16.3								0.0093															
17		0.0186	0.0944			0.0785		0.0851		0.1902	0.0638		0.1848		0.0279		0.0120			0.0745			0.2354
17.2							0.0040											0.0013					

Figure 43 United States—Asian-American population (minimum allele frequency = 0.0066)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
17.3								0.0532															
18		0.0013	0.2367			0.0372		0.0106		0.0319	0.1144		0.0891		0.1955		0.0013	0.0306		0.0718			0.1822
18.2							0.0013								0.0013								
18.3								0.0199															
19			0.1862			0.0399		0.0027		0.0013	0.1582		0.0053		0.1383			0.0346		0.0505			0.1130
19.3			0.0013					0.0053															
20			0.1662			0.0359					0.0931				0.0386			0.0612		0.0346			0.0146
21			0.1011			0.0239					0.0399				0.0040			0.1130		0.0239			0.0027
21.1						0.0013																	
21.3															0.0027								
22			0.1051			0.0213					0.0492				0.0013			0.1862		0.0239			
23			0.0532			0.0013					0.2168							0.2420		0.0080			
23.2																		0.0080					
24			0.0133			0.0040					0.1622							0.1569		0.0027			
24.2																		0.0066					
25			0.0120			0.0040					0.0745							0.1064		0.0013			
26			0.0027								0.0160							0.0426					
27			0.0013						0.0013		0.0027							0.0066					
28									0.0532		0.0013							0.0027					
28.2									0.0080														
29									0.2620														
29.2									0.0013														
30									0.2553														
30.2									0.0160														
30.3									0.0066														
31									0.1144														
31.2									0.0798														
32									0.0279														
32.2									0.1157														
33									0.0027														

United States—Asian-American population (continued)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA	
33.2									0.0532															
34.2									0.0013															
35.2									0.0013															

United States—Asian-American population (*continued*)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
Sample Size	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337	337
2.2																			0.0015				
5																				0.0682	0.0015		
6																			0.0015	0.0030	0.2255	0.0015	
7				0.0015												0.0119			0.0074	0.1528	0.1780		
8	0.0030			0.1098	0.0148									0.0074		0.1588	0.0193		0.0089	0.0119	0.1098	0.5059	
9	0.0267			0.0742	0.1291							0.0059		0.0534		0.1662	0.0134		0.1855	0.0148	0.1706	0.1276	
9.3																					0.3056		
10	0.2849			0.0682	0.0415	0.0119	0.0015	0.0030		0.0030		0.1988		0.0534	0.0074	0.2745	0.1083		0.1350	0.1068	0.0089	0.0490	
11	0.3056	0.0059		0.2834	0.3190	0.0089		0.0623		0.1380		0.3412	0.0030	0.3442	0.3086	0.2166	0.0653		0.1602	0.1246		0.2789	
11.3												0.0504											
12	0.3131	0.0341		0.3071	0.3056	0.1528	0.0653	0.1602		0.0045		0.0401		0.3724	0.2567	0.1454	0.1558		0.2181	0.1780		0.0371	
12.1								0.0015															
12.2								0.0015															
12.3												0.0015											
13	0.0638	0.2982		0.1113	0.1706	0.1187	0.2760	0.0697		0.0074	0.0015	0.0326	0.0015	0.1484	0.0742	0.0237	0.3249		0.2062	0.0831			0.0015
13.2							0.0163																
14	0.0030	0.2997		0.0445	0.0178	0.1751	0.3501	0.0623		0.0371	0.0015	0.2849	0.1528	0.0178	0.0490	0.0015	0.1914		0.0579	0.0519			0.0861
14.1																			0.0015				
14.2							0.0223																
14.3								0.0030															
15		0.1914	0.0445		0.0015	0.1528	0.1662	0.1528		0.3650	0.0015	0.0430	0.2774	0.0030	0.0134	0.0015	0.0905		0.0134	0.0475			0.1246
15.2							0.0356																
15.3								0.0890															
16		0.1320	0.0341			0.1098	0.0549	0.0979		0.3680	0.0430	0.0015	0.2329		0.0045		0.0282	0.0015	0.0030	0.0519			0.2166
16.2							0.0030																
16.3								0.0490															
17		0.0356	0.0994			0.1098	0.0030	0.0445		0.0712	0.1869		0.1988		0.0653		0.0030	0.0015		0.0415			0.2760
17.1								0.0030															

Figure 44 United States—Caucasian population (minimum allele frequency = 0.0074)



Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
17.2							0.0015																
17.3			0.0208					0.1246															
18		0.0030	0.1632			0.0861	0.0015	0.0030		0.0059	0.0846		0.1187		0.0638			0.0119		0.0326			0.1840
18.3			0.0223					0.0579															
19			0.1202			0.0445					0.1454		0.0148		0.1039			0.0549		0.0119			0.0979
19.3			0.0059					0.0178															
20			0.0994			0.0134					0.1573				0.0341			0.1513		0.0148			0.0134
20.2																		0.0045					
20.3			0.0015																				
21			0.1365			0.0104					0.0282				0.0119			0.1810		0.0015			
21.2																		0.0030					
21.3															0.0045								
22			0.1068			0.0030					0.0178				0.0015			0.1899					
22.2																		0.0089					
23			0.0816			0.0030					0.0979							0.1469					
23.2																		0.0045					
23.3															0.0015								
24			0.0386								0.1009							0.1454		0.0015			
25			0.0208								0.1128							0.0697		0.0015			
26			0.0030						0.0030		0.0178							0.0178					
27			0.0015						0.0282		0.0030							0.0059					
28									0.1632									0.0015					
29									0.2389														
29.2									0.0015														
29.3									0.0015														
30									0.2389														
30.2									0.0267														
31									0.0682														
31.2									0.0890														
32									0.0237														

United States—Caucasian population (continued)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
32.2									0.0920														
33.2									0.0208														
34.2									0.0045														

United States—Caucasian population (*continued*)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
Sample Size	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382	382
2.2																			0.0223				
3.2																			0.0026				
5																			0.0079	0.0524			
6					0.0013															0.0013	0.2788	0.0079	
7	0.0092		0.0026											0.0537		0.0118			0.0052	0.1021	0.3207	0.0013	
8	0.0052	0.0013		0.0916	0.0183									0.0131		0.1139	0.0065		0.0406	0.0380	0.0903	0.4830	
9	0.0236	0.0013		0.1597	0.1047	0.0013		0.0013				0.0013	0.0013	0.0484		0.0798	0.0026		0.1767	0.0118	0.1230	0.0733	
9.3			0.0013																		0.1715		
10	0.2605	0.0013		0.0942	0.1558	0.0065	0.0039	0.0026		0.0052		0.3089		0.0471	0.0105	0.2579	0.0982		0.1832	0.0694	0.0144	0.0563	
10.3																0.0013							
11	0.2814	0.0026		0.2277	0.3207	0.0118	0.0144	0.0366		0.0746		0.3220		0.3822	0.1649	0.2906	0.0510		0.1296	0.0668		0.2631	0.0013
11.1	0.0013																						
11.2							0.0026																
11.3												0.0432				0.0013							
12	0.3691	0.0445		0.2866	0.2435	0.1086	0.0812	0.0916		0.0105		0.0419	0.0013	0.2958	0.1715	0.1963	0.1257		0.1819	0.1688	0.0013	0.1086	0.0026
12.2							0.0144																
12.3												0.0039											
13	0.0419	0.2592	0.0013	0.0929	0.1401	0.1178	0.1872	0.0681		0.0092		0.0170	0.0039	0.1492	0.1047	0.0419	0.3220	0.0013	0.1688	0.1008		0.0013	0.0013
13.2							0.0707																
14	0.0052	0.3547	0.0013	0.0432	0.0131	0.1558	0.3010	0.1099		0.0223		0.2264	0.0877	0.0105	0.1427	0.0052	0.2356		0.0589	0.0707		0.0013	0.0654
14.2						0.0013	0.0445																
14.3								0.0026															
15	0.0026	0.2291	0.0366	0.0013	0.0026	0.1257	0.1309	0.1636		0.4254		0.0327	0.3442		0.0445		0.1178		0.0183	0.0812			0.0969
15.2						0.0013	0.0733																
15.3								0.0275															
16		0.0812	0.0497			0.1257	0.0406	0.1571		0.3599	0.0366	0.0026	0.2723		0.0065		0.0340		0.0013	0.0602			0.3076
16.1								0.0026															
16.2							0.0249																

Figure 45 United States—Hispanic population (minimum allele frequency = 0.0065)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
16.3								0.0537															
17		0.0236	0.0733	0.0026		0.1636	0.0065	0.0707		0.0798	0.1780		0.1780		0.0537		0.0065			0.0576			0.2709
17.1			0.0026																				
17.2							0.0039																
17.3			0.0118					0.1558															
18			0.2003			0.0851		0.0065		0.0118	0.0681		0.1021		0.0995			0.0065		0.0340			0.1819
18.3			0.0209					0.0432															
19			0.1885			0.0380					0.1846		0.0092		0.0720			0.0694	0.0026	0.0183			0.0654
19.3			0.0105					0.0065							0.0039								
20			0.1688			0.0183				0.0013	0.1361				0.0432			0.0903		0.0301			0.0065
20.2																		0.0026					
20.3															0.0118								
21			0.0916			0.0236					0.0366				0.0105			0.1401		0.0209			
21.3															0.0432								
22			0.0668			0.0065					0.0628							0.1387		0.0105			
22.2																		0.0065					
22.3															0.0131								
23			0.0366			0.0052					0.1401							0.1243		0.0052		0.0039	
23.2																		0.0039					
23.3															0.0039								
24		0.0013	0.0183			0.0026					0.0838							0.1649					
24.2									0.0026														
25			0.0118			0.0013					0.0550							0.1401					
26			0.0026						0.0039		0.0157							0.0681					
27			0.0026						0.0144		0.0013							0.0301					
28									0.1139		0.0013							0.0079					
28.2									0.0013														
29									0.2107									0.0039					
30									0.2723									0.0013					
30.2									0.0144														

United States—Hispanic population (continued)

Allele	CSF1PO	D10S1248	D12S391	D13S317	D16S539	D18S51	D19S433	D1S1656	D21S11	D22S1045	D2S1338	D2S441	D3S1358	D5S818	D6S1043	D7S820	D8S1179	FGA	PentaD	PentaE	TH01	TPOX	vWA
31									0.0497														
31.2									0.1139														
32									0.0157														
32.2									0.1257														
33									0.0013														
33.2									0.0550														
35									0.0026														
36									0.0013														
38									0.0013														

United States—Hispanic population (continued)

**Table 15 Random Match Probability (RMP) values for Verifiler™ Plus kit markers**

Allele	Combined China	African-American	Caucasian	Asian-American	Hispanic
CSF1PO	0.1168	0.0792	0.1289	0.1221	0.1289
D10S1248	0.0938	0.0672	0.0931	0.1007	0.0931
D12S391	0.0427	0.0362	0.0198	0.0408	0.0198
D13S317	0.0643	0.1363	0.0735	0.0675	0.0735
D16S539	0.0799	0.0727	0.0981	0.0836	0.0981
D18S51	0.0343	0.0300	0.0289	0.0366	0.0289
D19S433	0.0526	0.0361	0.0895	0.0549	0.0895
D1S1656	0.0440	0.0315	0.0195	0.0490	0.0195
D21S11	0.0509	0.0419	0.0466	0.0508	0.0466
D22S1045	0.0883	0.0504	0.1367	0.0844	0.1367
D2S1338	0.0318	0.0198	0.0279	0.0321	0.0279
D2S441	0.0858	0.0996	0.0976	0.0897	0.0976

Table 15 Random Match Probability (RMP) values for Verifiler Plus kit markers (continued)

Allele	Combined China	African-American	Caucasian	Asian-American	Hispanic
D3S1358	0.1262	0.1033	0.0757	0.1222	0.0757
D5S818	0.0870	0.0952	0.1290	0.0907	0.1290
D6S1043	0.0287	0.0248	0.0584	0.0313	0.0584
D7S820	0.0799	0.0747	0.0678	0.0905	0.0678
D8S1179	0.0449	0.0700	0.0604	0.0462	0.0604
FGA	0.0345	0.0301	0.0370	0.0393	0.0370
Penta D	0.0594	0.0237	0.0530	0.0577	0.0530
Penta E	0.0115	0.0188	0.0199	0.0121	0.0199
TH01	0.1490	0.0885	0.0811	0.1641	0.0811
TPOX	0.2129	0.0860	0.1785	0.2141	0.1785
vWA	0.0674	0.0602	0.0624	0.0641	0.0624
<b>Cumulative</b>	<b>1.34 x 10<sup>-28</sup></b>	<b>2.13 x 10<sup>-30</sup></b>	<b>1.24 x 10<sup>-28</sup></b>	<b>3.09 x 10<sup>-28</sup></b>	<b>1.24 x 10<sup>-28</sup></b>

## Probability of paternity exclusion

### Probability of paternity exclusion definition

The probability of paternity exclusion (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 STR loci in the kit (Chakraborty, Stivers, and Zhong, 1996).

### Probability of paternity exclusion observation

If your laboratory requires the probability of paternity exclusion calculation, contact HID Support.



# Troubleshooting

Observation	Possible cause	Recommended action
Faint or no signal from DNA Control 007 and the test samples at all loci, including the IQC markers	The incorrect volume of master mix or primer set was used.	Use the correct volume of master mix or primer set.
	The DNA polymerase was not activated.	Repeat the amplification with an initial hold for 1 minute at 95°C.
	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.
	The thermal cycler malfunctioned.	See the thermal cycler user guide and check the instrument calibration.
	Incorrect thermal cycler conditions were used.	Use the correct thermal cycler conditions.
	The tubes or plate were not seated tightly in the thermal cycler during amplification.	Push the tubes or plate firmly into the block after the first cycle.
	The wrong PCR reaction tubes or plate were used.	Use MicroAmp™ Reaction Tubes with Caps or a MicroAmp™ Optical 96-Well Reaction Plate.
	Insufficient PCR product was electrokinetically injected.	Use the correct settings for the capillary electrophoresis instrument.
Degraded formamide was used.	Ensure that the formamide is correctly stored. Do not thaw and refreeze the formamide multiple times. Try using Hi-Di™ Formamide.	
Positive signal from DNA Control 007 but partial or no signal from the 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 17.5 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 the test samples when IQC peaks are present and balanced <i>(continued)</i>	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, re-amplify 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 nuclease-free water or low-TE buffer (with 0.1 mM EDTA).
Positive signal from DNA Control 007 but partial or no signal from the test samples when IQC peaks are present and unbalanced	The test sample contains a high concentration of PCR inhibitor (for example, heme compounds, certain dyes).	Quantify the DNA, then use the minimum required volume of test sample DNA. Wash the sample in a Centricon™-100 centrifugal filter unit.
	The test sample was diluted in the wrong buffer (for example, a TE buffer with an incorrect EDTA concentration).	Redilute DNA using nuclease-free water or low-TE buffer (with 0.1 mM EDTA).
Positive signal from DNA Control 007 and elevated signal from the test samples when IQC peaks are present and unbalanced	The quantity of the test sample DNA is more than the 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 you apply the provided locus-specific stutter filters. See “Extra peaks: Stutter” on page 76.
	Stutter filters apply to sample input amounts $\geq 250$ pg. Samples well below the recommended input amount may exhibit stutter values that exceed the filters due to the stochastic effects of the PCR.	Increase the sample input to $>250$ pg, if possible. 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.

Observation	Possible cause	Recommended action
More than two alleles present at a locus (continued)	Incomplete 3' A base addition (n-1 bp position) occurred.	In the PCR, include the final extension step: 5 minutes 60°C.
		Remove the amplified plate from storage (thaw if needed) and place on the thermal cycler for 15 minutes at 60°C.
		Check the quantity of the original sample DNA to confirm that the input is <750 pg per reaction. If needed, adjust the input amount during re-amplification.
		If the total amount of DNA in the reaction is >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.	Ensure that you are using the recommended number of PCR cycles. Decrease the input DNA amount and repeat the PCR amplification, or interpret the off-scale data according to your laboratory procedure.
		Ensure 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 29 PCR cycles.	
The double-stranded DNA was not completely denatured.	Use the recommended amount of Hi-Di™ Formamide and heat the sample plate for 3 minutes at 95°C.	
Poor peak height balance	Incorrect thermal cycler conditions were used.	Use the correct thermal cycler conditions.
Some, but not all, loci are visible on test sample electropherograms	The DNA quantity was too low, leading to stochastic effects.	Load more DNA sample or concentrate the DNA.
STR profiles contain many off-scale alleles	DNA quantification was not performed or was not accurate.	Ensure that DNA quantification is accurate.



# Materials required but not supplied

- Sample preparation (PCR or CE) required materials ..... 123
- Thermal cycler required materials ..... 123
- Capillary electrophoresis instrument required materials ..... 124
- Analysis software required materials ..... 126
- Miscellaneous required materials ..... 126

Unless otherwise indicated, all materials are available through [thermofisher.com](http://thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](http://fisherscientific.com) or another major laboratory supplier.

Catalog numbers that appear as links open the web pages for those products.

## Sample preparation (PCR or CE) required materials

Item	Amount	Cat. No.
GeneScan™ 600 LIZ™ Size Standard v2.0 <b>IMPORTANT!</b> Do not use GeneScan™ 350 ROX™, GeneScan™ 500 ROX™, or GeneScan™ 500 LIZ™ Size Standards with this kit.	2 × 200 µL	<a href="#">4408399</a>
TE Buffer [low-TE buffer; 10 mM Tris-HCl (pH 8.0) and 0.1 mM EDTA]	100 mL	<a href="#">12090015</a> or see "(Optional) Prepare low-TE buffer" on page 20
Nuclease-Free Water (not DEPC-Treated)	10 × 50 mL	<a href="#">AM9937</a>
Hi-Di™ Formamide	25 mL	<a href="#">4311320</a>

## Thermal cycler required materials

### ProFlex™ PCR System

Item	Source
ProFlex™ 96-well PCR System	<a href="#">4484075</a>
ProFlex™ 2 × 96-well PCR System	<a href="#">4484076</a>
ProFlex™ 3 × 32-Well PCR System	<a href="#">4484073</a>



## VeritiPro™ Thermal Cycler

Item	Source
HID VeritiPro™ Thermal Cycler, 96-well	<a href="#">A52127</a>

## Capillary electrophoresis instrument required materials

### SeqStudio™ Flex Series Genetic Analyzer for Human Identification

Item	Source
SeqStudio™ 8 Flex Genetic Analyzer for Human Identification with SeqStudio™ Flex Series Instrument Software v1.1.1	<a href="#">A56532</a>
SeqStudio™ 24 Flex Genetic Analyzer for Human Identification with SeqStudio™ Flex Series Instrument Software v1.1.1	<a href="#">A56534</a>
Anode Buffer Container 3500/Flex Series	<a href="#">4393927</a>
Cathode Buffer Container 3500/Flex Series	<a href="#">4408256</a>
Septa Cathode Buffer Container 3500/Flex Series	<a href="#">4410715</a>
Capillary array 36-cm SeqStudio™ 8 Flex	<a href="#">A49104</a>
Capillary array 36-cm SeqStudio™ 24 Flex	<a href="#">A49105</a>
96-Well Standard Retainer & Base Set SeqStudio™ Flex Series	<a href="#">A49316</a>
8-Tube Standard Retainer & Base Set SeqStudio™ Flex Series	<a href="#">A49296</a>
8-Strip Septa 3500/Flex Series	<a href="#">4410701</a>
Septa for 96-Well Plates, for 3500/SeqStudio™ Flex	<a href="#">4412614</a>
DS-37 Matrix Standard Kit (Dye set J6-T, 6-dye)	<a href="#">A31234</a>
POP-4™ (960) Performance Optimized Polymer	<a href="#">4393710</a>
POP-4™ (384) Performance Optimized Polymer	<a href="#">4393715</a>
Conditioning Reagent Kit 3500/Flex Series	<a href="#">4393718</a>

## SeqStudio™ Genetic Analyzer for HID

Item	Source
SeqStudio™ Genetic Analyzer for HID with SeqStudio™ Data Collection Software v1.2.1, v1.2.4, or v1.2.5	<a href="#">A46227</a>
(Optional) SAE Administrator Console v2.0 or v2.1	<a href="#">A46170</a> or <a href="#">A53717</a>
(Optional) SeqStudio™ Plate Manager v1.2 or v1.3	Available on <a href="https://apps.thermofisher.com">apps.thermofisher.com</a> or for download at <a href="https://thermofisher.com">thermofisher.com</a>
SeqStudio™ Genetic Analyzer Cartridge v2	<a href="#">A41331</a>
SeqStudio™ Genetic Analyzer Cathode Buffer Container	<a href="#">A33401</a>
Reservoir Septa (for Cathode Buffer Container)	<a href="#">A35640</a>
SeqStudio™ Integrated Capillary Protector	<a href="#">A31923</a>
Septa for SeqStudio™ Genetic Analyzer, 96 well	<a href="#">A35641</a>
Septa for SeqStudio™ Genetic Analyzer, 8 strip	<a href="#">A35643</a>
DS-37 Matrix Standard Kit (Dye set J6-T, 6-dye)	<a href="#">A31234</a>

## 3500 Series Genetic Analyzer for Human Identification

Item	Source
3500 Genetic Analyzer for Human Identification with 3500 Series Data Collection Software 4	<a href="#">4406017</a>
3500xL Genetic Analyzer for Human Identification with 3500 Series Data Collection Software 4	<a href="#">4406016</a>
(Software upgrade package) 3500 Series HID Data Collection Software v4.0.1	<a href="#">A46085</a>
Anode Buffer Container 3500/Flex Series	<a href="#">4393927</a>
Cathode Buffer Container 3500/Flex Series	<a href="#">4408256</a>
Septa Cathode Buffer Container 3500/Flex Series	<a href="#">4410715</a>
3500 Genetic Analyzer 8-Capillary Array, 36 cm	<a href="#">4404683</a>
3500xL Genetic Analyzer 24-Capillary Array, 36 cm	<a href="#">4404687</a>
Retainer & Base Set (Standard) for 3500/3500xL Genetic Analyzer, 96 well	<a href="#">4410228</a>
Retainer & Base Set (Standard) for 3500/3500xL Genetic Analyzer, 8 tube	<a href="#">4410231</a>
8-Strip Septa 3500/Flex Series	<a href="#">4410701</a>
Septa for 96-Well Plates, for 3500/SeqStudio™ Flex	<a href="#">4412614</a>
DS-37 Matrix Standard Kit (Dye set J6-T, 6-dye)	<a href="#">A31234</a>
POP-4™ (960) Performance Optimized Polymer	<a href="#">4393710</a>

(continued)

Item	Source
POP-4™ (384) Performance Optimized Polymer	4393715
Conditioning Reagent Kit 3500/Flex Series	4393718

## Analysis software required materials

### GeneMapper™ ID-X Software

Item	Source
GeneMapper™ ID-X Software v1.7.2 patch <sup>[1]</sup>	Thermo Fisher Scientific <sup>[2]</sup>
GeneMapper™ ID-X Software v1.7 Full Installation	A71700
GeneMapper™ ID-X Software v1.7 Client Installation	A71701
GeneMapper™ ID-X Software v1.6 Full Installation	A39975
GeneMapper™ ID-X Software v1.6 Client Installation	A39976
GeneMapper™ ID-X Software v1.5 Full Installation	A27884
GeneMapper™ ID-X Software v1.5 Client Installation	A27886

<sup>[1]</sup> The patch addresses known issues and provides new user functionality since the v1.7 release.

<sup>[2]</sup> Available for free download at [www.thermofisher.com/GMIDXsoftware](http://www.thermofisher.com/GMIDXsoftware).

---

**Note:** For a list of GeneMapper™ ID-X Software versions that are compatible with your kit and capillary electrophoresis instrument, see “Instruments and software compatibility” on page 15.

---

## 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-Tube Strip, 0.2 mL	4316567
MicroAmp™ Optical 8-Cap Strips	4323032

(continued)

Item	Source
MicroAmp™ 96-Well Tray/Retainer Set (Adapter for 8-Tube Strip)	403081
MicroAmp™ 96-Well Base	N8010531
MicroAmp™ Clear Adhesive Film	4306311
MicroAmp™ Optical Adhesive Film	4311971
MicroAmp™ Optical 96-Well Reaction Plate	N8010560
MicroAmp™ Optical 96-Well Reaction Plate with Barcode	4326659

## Laboratory supplies

Item	Source
<b>Various procedures</b>	
Aerosol resistant pipette tips	MLS <sup>[1]</sup>
Microcentrifuge tubes	MLS
Pipettors	MLS
Tape, labeling	MLS
Tube, 50-mL Falcon™	MLS
Tube decapper, autoclavable	MLS
Deionized water, PCR grade	MLS
Vortex	MLS
(Optional) Tabletop centrifuge with 96-Well Plate Adapters	MLS
(Optional) Handheld Barcode Scanner	4488442

<sup>[1]</sup> Major laboratory supplier



# PCR work areas

■ Work area setup and lab design .....	128
■ PCR setup work area materials .....	128
■ Amplified DNA work area .....	129

## 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
- Pipettes





- Tube decapper that can be autoclaved
- 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, visit [thermofisher.com/support](https://www.thermofisher.com/support).

## 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 sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer 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 needed) 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.



**AVERTISSEMENT ! PRÉCAUTIONS GÉNÉRALES EN CAS DE MANIPULATION DE PRODUITS CHIMIQUES.** Pour minimiser les risques, veiller à ce que le personnel du laboratoire lise attentivement et mette en œuvre les consignes de sécurité générales relatives à l'utilisation et au stockage des produits chimiques et à la gestion des déchets qui en découlent, décrites ci-dessous. Consulter également la FDS appropriée pour connaître les précautions et instructions particulières à respecter :

- Lire et comprendre les fiches de données de sécurité (FDS) fournies par le fabricant avant de stocker, de manipuler ou d'utiliser les matériaux dangereux ou les produits chimiques. Pour obtenir les FDS, se reporter à la section « Documentation et support » du présent document.
- Limiter les contacts avec les produits chimiques. Porter des équipements de protection appropriés lors de la manipulation des produits chimiques (par exemple : lunettes de sûreté, gants ou vêtements de protection).
- Limiter l'inhalation des produits chimiques. Ne pas laisser les récipients de produits chimiques ouverts. Ils ne doivent être utilisés qu'avec une ventilation adéquate (par exemple, sorbonne).
- Vérifier régulièrement l'absence de fuite ou d'écoulement des produits chimiques. En cas de fuite ou d'écoulement d'un produit, respecter les directives de nettoyage du fabricant recommandées dans la FDS.
- Manipuler les déchets chimiques dans une sorbonne.

- Veiller à utiliser des récipients à déchets primaire et secondaire. (Le récipient primaire contient les déchets immédiats, le récipient secondaire contient les fuites et les écoulements du récipient primaire. Les deux récipients doivent être compatibles avec les matériaux mis au rebut et conformes aux exigences locales, nationales et communautaires en matière de confinement des récipients.)
- Une fois le récipient à déchets vidé, il doit être refermé hermétiquement avec le couvercle fourni.
- Caractériser (par une analyse si nécessaire) les déchets générés par les applications, les réactifs et les substrats particuliers utilisés dans le laboratoire.
- Vérifier que les déchets sont convenablement stockés, transférés, transportés et éliminés en respectant toutes les réglementations locales, nationales et/ou communautaires en vigueur.
- **IMPORTANT !** Les matériaux représentant un danger biologique ou radioactif exigent parfois une manipulation spéciale, et des limitations peuvent s'appliquer à leur élimination.

## Biological hazard safety



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



**WARNING! BIOHAZARD.** Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. 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)*, 6th Edition, HHS Publication No. (CDC) 300859, Revised June 2020  
[cdc.gov/labs/bmbi](https://www.cdc.gov/labs/bmbi)
- Laboratory biosafety manual, fourth edition. Geneva: World Health Organization; 2020 (Laboratory biosafety manual, fourth edition and associated monographs)  
[who.int/publications/i/item/9789240011311](https://www.who.int/publications/i/item/9789240011311)



## Related documentation

Table 16 STR and quantification kits

Document title	Pub. No.
<b>STR kits</b>	
<i>Verifiler™ Plus PCR Amplification Kit—PCR Amplification and CE Quick Reference</i>	MAN0017495
<i>Verifiler™ Plus PCR Amplification Kit—PCR Setup Quick Reference</i>	MAN0017494
<i>Technical Note: Direct Amplification of Reference Samples Using the Verifiler™ Plus PCR Amplification Kit</i>	Go to <a href="https://www.thermofisher.com">thermofisher.com</a> , then search for the technical note by title, or contact your local Human Identification representative.
<i>Technical Note: Handling STR Kits and Ladder Decontamination</i>	
<b>Quantification kits</b>	
<i>Quantifiler™ HP and Quantifiler™ Trio DNA Quantification Kits User Guide</i>	4485355

Table 17 Thermal cyclers

Document title	Pub. No.
<b>ProFlex™ PCR System</b>	
<i>ProFlex™ PCR System User Guide</i>	<a href="#">MAN0007697</a>
<i>ProFlex™ PCR System Kit Validation User Bulletin</i>	100031595 <sup>[1]</sup>
<b>VeritiPro™ Thermal Cycler</b>	
<i>VeritiPro™ Thermal Cycler User Guide</i>	<a href="#">MAN0019157</a>
<i>HID VeritiPro™ Thermal Cycler, 96-well, User Bulletin—Applied Biosystems™ PCR Amplification Kit Validation</i>	<a href="#">MAN0025561</a>

<sup>[1]</sup> Archived document. To access, go to [https://assets.thermofisher.com/TFS-Assets/LSG/manuals/100031595\\_ProFlexKit\\_Validation\\_UB.pdf](https://assets.thermofisher.com/TFS-Assets/LSG/manuals/100031595_ProFlexKit_Validation_UB.pdf)



**Table 18 Capillary electrophoresis instruments**

Document title	Pub. No.
<b>SeqStudio™ Flex Series Genetic Analyzer for Human Identification</b>	
<i>SeqStudio™ Flex Series Genetic Analyzer with Instrument Software v1.1.1 User Guide</i>	100104689
<i>SeqStudio™ Flex Series Genetic Analyzer for HID Validation User Bulletin</i>	MAN0028463
<i>SeqStudio™ Flex Series Instrument Software v1.1.1 User Bulletin</i>	MAN0029757
<b>SeqStudio™ Genetic Analyzer for HID</b>	
<i>SeqStudio™ Genetic Analyzer Instrument and Software User Guide (v1.2 and later)</i>	MAN0018646
<i>SeqStudio™ Genetic Analyzer for HID User Bulletin—New Software Features and Verification/Validation Studies (v1.2 and later)</i>	MAN1001221
<b>3500 Series Genetic Analyzer for Human Identification</b>	
<i>3500 Series Data Collection Software 4 User Bulletin: New Features and Developmental Validation</i>	100075298
<i>3500/3500xL Genetic Analyzer with 3500 Series Data Collection Software v3.3 User Guide</i>	100079380

**Table 19 Analysis software**

Document title	Pub. No.
<b>GeneMapper™ ID-X Software all versions</b>	
<i>GeneMapper™ ID-X Software Bin Overlap User Bulletin</i>	100029546
<i>Technical Note: Customizing GeneMapper™ ID-X Software Panel and Bin Sets to Include or Exclude Internal Quality Control Markers in the Quality Value System (v1.5.2 and later)</i>	Thermo Fisher Scientific <sup>[1]</sup>
<i>Technical Note: Compendium of GeneMapper™ ID-X Software version changes from version 1.0.1 through version 1.7.2</i>	
<b>GeneMapper™ ID-X Software v1.7</b>	
<i>GeneMapper™ ID-X Software v1.7 Administration Guide</i>	MAN0029245
<i>GeneMapper™ ID-X Software v1.7 Installation Guide</i>	MAN0029246
<i>GeneMapper™ ID-X Software v1.7 New Features and Software Verification and Validation User Bulletin</i>	MAN0029209
<b>GeneMapper™ ID-X Software v1.6</b>	
<i>GeneMapper™ ID-X Software v1.6 New Features and Software Verification User Bulletin</i>	100073905
<b>GeneMapper™ ID-X Software v1.5</b>	
<i>GeneMapper™ ID-X Software v1.5 New Features and Verification User Bulletin</i>	100031708
<i>GeneMapper™ ID-X Software v1.5 Getting Started Guide— Basic Features</i>	100031701

**Table 19 Analysis software** (continued)

Document title	Pub. No.
<i>GeneMapper™ ID-X Software v1.5 Quick Reference— Basic Features</i>	100031702
<i>GeneMapper™ ID-X Software v1.5 Getting Started Guide— Mixture Analysis Tool</i>	100031704
<i>GeneMapper™ ID-X Software v1.5 Quick Reference— Mixture Analysis Tool</i>	100031705
<i>GeneMapper™ ID-X Software v1.5 Installation Guide</i>	100031706
<i>GeneMapper™ ID-X Software v1.5 Administrator Guide</i>	100031703
<i>GeneMapper™ ID-X Software v1.5 Reference Guide</i>	100031707

<sup>[1]</sup> Go to [thermofisher.com](http://thermofisher.com), then search for the technical note by title, or contact your local Human Identification representative.

## Customer and technical support

For support, use one of the contact methods listed below, depending on your location.

Location	Contact method
In North America	Send an email to: <a href="mailto:HIDTechSupport@thermofisher.com">HIDTechSupport@thermofisher.com</a>
	Call 888-821-4443; select option 2, say "Application Support", then say "HID" or "Human Identification".
Outside North America	Contact your local support office.

For the latest services and support information for all locations, go to [thermofisher.com/support](http://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 its affiliates 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](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://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/](http://www.springerlink.com/content/p713q3w5440674u3/)
- Burri, H., Sulzer, A., Voegeli, P. and Kratzer, A. 2015. Accounting for linkage between the STR loci D5S818/CSF1PO and vWA/D12S391 in kinship analyses: Impact on likelihood ratio values. *Forensic Sci. Int.* 5:E387–E389. (<http://dx.doi.org/10.1016/j.fsigss.2015.09.153>)
- Butler, J.M. 2005. *Forensic DNA Typing*. Burlington, MA:Elsevier Academic Press.
- Butler, J.M. 2010. *Fundamentals of 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. Arlequin suite v. 3.5. *Mol. Ecol. Res.* 10:564–567.



- Frank, W., Llewellyn, B., Fish, P., et al. 2001. Validation of the AmpFISTR 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., Matisse, 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 AmpFISTR NGM SElect PCR Amplification Kit for Han population in Fujian Province, China. *Int J. Legal Med.* 127:345–346.
- Liu, Q.-L., Luo, H., Zhao, H., Huang, X.-L., Cheng, J.-D. and Lu, D.-J. 2013. Recombination analysis of autosomal short tandem repeats in Chinese Han families. *Electrophoresis* 35(6):883–887. (<https://doi.org/10.1002/elps.201300417>)
- 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]<sub>n</sub> 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 reimplement of the Genepop software for Windows and Linux. *Molecular Ecology Resources* 8:103–106.
- Sensabaugh, G.F. 1982. Biochemical markers of individuality. In: Saferstein, R., ed. *Forensic Science Handbook*. Prentice-Hall, Inc., New York, pp. 338–415.
- Sharma, V. and Litt, M. 1992. Tetranucleotide repeat polymorphism at the D21S11 locus. *Hum Mol. Genet.* 1:67.
- 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 (SWGDM). 2016. Validation Guidelines for DNA Analysis Methods. Available at [http://media.wix.com/ugd/4344b0\\_813b241e8944497e99b9c45b163b76bd.pdf](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 AmpFISTR 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.



