

AmpF \mathcal{L} STR® SEfiler Plus™ PCR Amplification Kit

Publication Number 4382731 Rev. D Revision Date August 2012

IMPORTANT! Reagents used for amplification must be isolated from any source of contaminating DNA, especially from previously amplified AmpF \mathcal{L} STR® PCR products and the AmpF \mathcal{L} STR® Allelic Ladders.

Thaw reagents completely before first use. To avoid freeze/thaw cycles, store components at 2 to 8°C after initial use.

The AmpF \mathcal{L} STR® SEfiler Plus™ Primer Set and Allelic Ladder contain primers that are labeled with light-sensitive dyes. Protect the primers and allelic ladder from light; do not leave them exposed on the laboratory bench.

To obtain optimal amplification and typing results, follow the procedures in the *AmpF \mathcal{L} STR® SEfiler Plus™ PCR Amplification Kit User Guide* (Pub. no. 4385739).

Table 1 AmpF \mathcal{L} STR® SEfiler Plus™ PCR Amplification Kit 200 reactions (Part no. 4382699)

Part no.	Part	Storage Conditions
4382711	AmpF \mathcal{L} STR® SEfiler Plus™ Master Mix, 2 tubes, 1.1 mL/tube Contains MgCl ₂ , dATP, dGTP, dCTP, and dTTP, bovine serum albumin, enzyme, and 0.05% sodium azide in buffer and salt.	-15 to -25°C upon receipt. 2 to 8°C after initial use.
4307984	AmpF \mathcal{L} STR® Control DNA 007, 1 tube, 0.3 mL Contains 0.10 ng/ μ L human male genomic DNA in 0.05% sodium azide and buffer. The profile of this male DNA is D2S1338 20, 23; D3S1358 15, 16; D8S1179 12, 13; D16S539 9, 10; D18S51 12,15; D19S433 14, 15; D21S11 28, 31; FGA 24, 26; SE-33 17, 25.2; TH01 7, 9.3; vWA 14, 16; Amelogenin X, Y.	
4382708	AmpF \mathcal{L} STR® SEfiler Plus™ Primer Set, 1 tube, 0.5 mL Contains locus-specific 6-FAM™, VIC®, NED™, and PET® dye-labeled and unlabeled primers in buffer. The primers amplify the STR loci D2S1338, D3S1358, D8S1179, D16S539, D18S51, D19S433, D21S11, FGA, SE-33, TH01, vWA, and the gender marker Amelogenin.	-15 to -25°C upon receipt. 2 to 8°C after initial use. Store protected from light.
4382703	AmpF \mathcal{L} STR® SEfiler Plus™ Allelic Ladder, 1 tube, 0.05 mL Contains the following amplified alleles: <ul style="list-style-type: none"> 6-FAM™ dye (blue): D3S1358 12-19; vWA 11-24; D16S539 5, 8-15; D2S1338 15-28. VIC® dye (green): Amelogenin X, Y; D8S1179 8-19; SE-33 4.2, 6.3, 8-37, 21.1, 21.2, 22.2, 23.2, 24.2, 25.2, 26.2, 27.2, 28.2, 29.2, 30.2, 31.2, 32.2, 33.2, 34.2, 35.2. NED™ dye (yellow): D19S433 9-17,12.2, 13.2, 14.2, 15.2, 16.2, 17.2; TH01 4-11, 9.3, 13.3; FGA 17-26, 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, 50.2, 51.2. PET® dye (red): D21S11 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, 35.2, 36-38; D18S51 7, 9, 10, 10.2, 11-13, 13.2, 14, 14.2, 15-27. 	-15 to -25°C upon receipt. 2 to 8°C after initial use. Store protected from light. IMPORTANT! The allelic ladder contains PCR products. Do NOT amplify. To avoid contamination, store the allelic ladder separate from the other kit components and unamplified DNA.

Note: For safety and biohazard guidelines, refer to the “Safety” section in the *AmpF \mathcal{L} STR® SEfiler Plus™ PCR Amplification Kit User Guide* (Pub. no. 4385739). For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Table 2 Materials required but not included in the kit



Part no.	Part
4366589	GeneScan™ 600 LIZ® Size Standard v2.0, 2 tubes, 200 µL/tube
4385739	AmpFSTR® SEfiler Plus™ PCR Amplification Kit User Guide

Performance Characteristics

The AmpFSTR® Kits are developed and manufactured by Life Technologies Corporation. We test each lot of kits to ensure that the kits perform according to specifications and are free from interfering contaminants.

Each AmpFSTR® SEfiler Plus™ PCR Amplification Kit contains reagents optimized to amplify and type 0.5 ng of AmpFSTR® Control DNA 007.

The recommended range of input sample DNA is approximately 0.5 to 0.75 ng of human DNA. However, we have successfully typed samples containing less than 0.5 ng of human DNA.

Follow the protocol in the *AmpFSTR® SEfiler Plus™ PCR Amplification Kit User Guide* to collect and analyze data on the instruments and software supported for use with this kit.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

For support visit www.lifetechnologies.com/support

www.lifetechnologies.com

