

GlobalFiler™ PCR Amplification Kit (1,000 Reactions)

Catalog Numbers 4482815, 4476135

Pub. No. 4477595 Rev. D

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *GlobalFiler™ PCR Amplification Kit User Guide* (Pub. No. 4477604). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Important guidelines

- Thaw reagents completely before first use. Store components at 2–8°C after first use to avoid freeze/thaw cycles.
- Keep reagents away from any source of contaminating DNA, especially from previously amplified PCR products and the allelic ladder.
- The 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 results, follow the procedures in the *GlobalFiler™ PCR Amplification Kit User Guide* (Pub. No. 4477604).

Contents and storage

Contents	Cat. No. 4482815 1,000 reactions	Storage
GlobalFiler™ Master Mix Contains MgCl ₂ , dATP, dGTP, dCTP, and dTTP, bovine serum albumin, enzyme, and 0.05% sodium azide in buffer and salt.	1 bottle, 7.5 mL	–25°C to –15°C on receipt. 2°C to 8°C after first use for up to 6 months or up to the expiration date stated on the kit (whichever comes first).
GlobalFiler™ Primer Set Contains locus-specific 6-FAM™, VIC™, NED™, TAZ™, and SID™ dye-labeled and unlabeled primers in buffer. The primers amplify the STR loci D3S1358, vWA, D16S539, CSF1PO, TPOX, D8S1179, D21S11, D18S51, DYS391, D2S441, D19S433, TH01, FGA, D22S1045, D5S818, D13S317, D7S820, SE33, D10S1248, D1S1656, D12S391, D2S1338, and the sex-determining markers, Y indel and Amelogenin.	1 bottle, 2.5 mL	–25°C to –15°C on receipt. 2°C to 8°C after first use for up to 6 months or up to the expiration date stated on the kit (whichever comes first). Store protected from light.

Contents	Cat. No. 4482815 1,000 reactions	Storage
GlobalFiler™ Allelic Ladder Contains the following amplified alleles: <ul style="list-style-type: none"> • 6-FAM™ dye (blue): D3S1358 9–20; vWA 11–24; D16S539 5, 8–15; CSF1PO 6–15; TPOX 5–15. • VIC™ dye (green): Y indel 1, 2; Amelogenin X, Y; D8S1179 5–19; 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, 35.2, 36–38; D18S51 7, 9, 10, 10.2, 11–13, 13.2, 14, 14.2, 15–27; DYS391 7–13. • NED™ dye (yellow): D2S441 8–11, 11.3, 12–17; D19S433 6–12, 12.2, 13, 13.2, 14, 14.2, 15, 15.2, 16, 16.2, 17, 17.2, 18.2, 19.2; TH01 4–9, 9.3, 10, 11, 13.3; FGA 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, 50.2, 51.2. • TAZ™ dye (red): D22S1045 8–19; D5S818 7–18; D13S317 5–16; D7S820 6–15; SE33 4.2, 6.3, 8, 9, 11–20, 20.2, 21, 21.2, 22.2, 23.2, 24.2, 25.2, 26.2, 27.2, 28.2, 29.2, 30.2, 31.2, 32.2, 33.2, 34.2, 35, 35.2, 36, 37. • SID™ dye (purple): D10S1248 8–19; D1S1656 9–14, 14.3, 15, 15.3, 16, 16.3, 17, 17.3, 18.3, 19.3, 20.3; D12S391 14–19, 19.3, 20–27; D2S1338 11–28. 	1 tube, 0.15 mL	–25°C to –15°C on receipt. 2°C to 8°C after first use or up to the expiration date stated on the kit. 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.
DNA Control 007 Contains 0.1 ng/μL human male genomic DNA in 0.05% sodium azide and buffer. The profile of this male DNA is D3S1358 15, 16; vWA 14, 16, D16S539 9, 10; CSF1PO 11, 12; TPOX 8; Y indel 2; Amelogenin X, Y; D8S1179 12, 13; D21S11 28, 31; D18S51 12, 15; DYS391 11; D2S441 14, 15; D19S433 14, 15; TH01 7, 9.3; FGA 24, 26; D22S1045 11, 16; D5S818 11; D13S317 11 D7S820 7, 12; SE33 17, 25.2; D10S1248 12, 15; D1S1656 13, 16; D12S391 18, 19; D2S1338 20, 23.	1 tube, 0.6 mL	–25°C to –15°C on receipt. 2°C to 8°C after first use or up to the expiration date stated on the kit.

Required materials not supplied

Unless otherwise indicated, all materials are available through thermofisher.com.

Item	Cat. No.
GeneScan™ 600 LIZ™ Size Standard v2.0, 2 tubes, 200 μL each IMPORTANT! Do not use GeneScan™ 350 ROX™, GeneScan™ 500 ROX™, or GeneScan™ 500 LIZ™ Size Standards with this kit.	4408399
Low-TE buffer (10 mM Tris, 0.1 mM EDTA, pH 8.0)	Teknova T0223

Performance characteristics

Each GlobalFiler™ PCR Amplification Kit is developed and manufactured by Thermo Fisher Scientific. We test each lot of kits to ensure that the kits perform according to specifications.

Each GlobalFiler™ PCR Amplification Kit contains reagents that are optimized to amplify and type 1 ng of DNA Control 007.

The recommended range of input sample DNA is approximately 0.5 ng to 1.5 ng of human DNA. However, we have successfully typed samples containing less than 0.5 ng of human DNA.

Follow the protocol in the *GlobalFiler™ PCR Amplification Kit User Guide* to collect and analyze data on the instruments and software that is 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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



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

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

Revision history: Pub. No. 4477595

Revision	Date	Description
D	19 October 2018	Updated branding and trademarks, no technical changes.
C	3 April 2015	Removed Master Mix Additive.
B	30 April 2014	Change storage condition from 2 months to 6 months. Update Globalfiler to a registered trademark.
A	3 June 2013	Initial publication.

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

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.