

Quantifiler™ Trio Automated DNA Quantification Kit

Catalog Number A58787

Pub. No. MAN1000067 Rev. A00

Note: For safety and biohazard guidelines, see the “Safety” appendix in the following product documentation: *Quantifiler™ Trio Automated DNA Quantification Kit User Guide* (Pub. No. MAN1000066). 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.
- The primer mix contains primers that are labeled with light-sensitive dyes. The reaction mix contains a light-sensitive dye. Protect the primer mix and the reaction mix from light; do not leave them exposed on the laboratory bench.
- To obtain optimal results, follow the procedures in the *Quantifiler™ Trio Automated DNA Quantification Kit User Guide* (Pub. No. MAN1000066).

Contents and storage

Table 1 Quantifiler™ Trio Automated DNA Quantification Kit (Cat. No. [A58787](#))

Item	Amount	Description	Storage ^[1]
Quantifiler™ THP PCR Reaction Mix	10 x 1 mL	Contains dNTPs, buffer, enzyme, MUSTANG PURPLE™ passive reference standard, and stabilizers.	–25°C to –15°C upon receipt. 2°C to 8°C after initial use. Protect from light. ^[2]
Quantifiler™ Trio Primer Mix	10 x 0.8 mL	Contains target-specific primers, ABY™, JUN™, VIC™, and FAM™ dye-labeled probes, and Internal PCR Control (IPC) template.	
Quantifiler™ THP DNA Dilution Buffer	6 x 1.8 mL	Contains genomic DNA standard dilution buffer.	–25°C to –15°C upon receipt. 2°C to 8°C after initial use.
Quantifiler™ THP DNA Standard	2 x 0.12 mL	Contains genomic DNA standard formulated at 100 ng/μL to generate standard curves.	
Quantifiler™ Automation Enhancer	1 x 0.2 mL	Contains defoaming agent.	15°C to 30°C upon receipt and after initial use.

^[1] See the expiration date on the package. Do not use expired product.

^[2] Excessive exposure to light can affect the fluorescent probes and the passive reference dye.

Performance characteristics

The Quantifiler™ Trio Automated DNA Quantification Kit is developed and manufactured by Thermo Fisher Scientific. We test each lot of kits to help ensure that the kits perform according to specifications.

Each kit contains reagents optimized to quantify DNA sample concentrations of 5 pg/μL to 100 ng/μL (depending on the range of DNA quantification standards used). Follow the protocols described in the *Quantifiler™ Trio Automated DNA Quantification Kit User Guide* (Pub. No. MAN1000066) and run the samples on the Applied Biosystems™ 7500 Real-Time PCR System and the Applied Biosystems™ QuantStudio™ 5 Real-Time PCR System.

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.

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

Revision history: Pub. No. MAN1000067 A00

Revision	Date	Description
A00	14 March 2024	New document for the Quantifiler™ Trio Automated DNA Quantification 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: 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.

©2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.