



Quantifiler™ Duo DNA Quantification Kit

Catalog Number 4387746

Pub. No. 4387913 Rev. F

Contents and storage

Contents	Quantity	Storage conditions
Quantifiler™ Duo PCR Reaction Mix Bottle of PCR Reaction mix containing dNTPs, buffer, AmpliTaq Gold™ DNA Polymerase, and ROX™ Passive Reference Standard and stabilizers.	5.0 mL	–15°C to –25°C upon receipt Store at 2°C to 8°C after first use
Quantifiler™ Duo DNA Standard Tube of genomic DNA Standard formulated at 200 ng/L to generate Standard Curves	0.12 mL	–15°C to –25°C upon receipt Store at 2°C to 8°C after first use
Quantifiler™ Duo Primer Mix Three tubes containing target-specific primers and FAM™, VIC™, and NED™ dye-labeled probes and Internal PCR Control (IPC) template.	1.4 mL/tube	–15°C to –25°C upon receipt Store at 2°C to 8°C after first use  CAUTION! The Quantifiler™ Duo Primer Mix contains probes that are labeled with light-sensitive dyes. Thaw completely at room temperature protected from light. Vortex 3 to 5 seconds before use.
Quantifiler™ Duo DNA Std. Dilution Buffer Two tubes containing DNA Standard dilution buffer.	1.8 mL/tube	–15°C to –25°C upon receipt Store at 2°C to 8°C after first use

 **CAUTION!** Reagents that are used for quantification should be isolated from contaminating sources of DNA, especially previously amplified products.

For safety and biohazard guidelines, see the Safety section in the *Quantifiler™ Duo DNA Quantification Kit User Guide* (Pub. No. 4391294). For the Quantifiler™ Duo PCR Reaction Mix, read the MSDS (4345050) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

To obtain optimal amplification and quantification results, it is critical that the protocols in the *Quantifiler™ Duo DNA Quantification Kit User Guide* (Pub. No. 4391294) are followed. The user guide is purchased separately.

Materials required but not supplied

Cat. No.	Material	Description
4391294	<i>Quantifiler™ Duo DNA Quantification Kit User Guide</i>	Included are all the necessary protocols for use with the Quantifiler™ Duo DNA Quantification Kit. Also included are protocols for setting up the Applied Biosystems™ 7500 Real-Time PCR System. To obtain optimal results, it is critical that the procedures detailed in the user guide are followed. Additional information on the results and their interpretation, a summary of experiments performed and results obtained, and troubleshooting information are also provided in the user guide.

Performance characteristics

The Quantifiler™ Duo DNA Quantification Kit is developed and manufactured by Thermo Fisher Scientific. Each lot of the kit is carefully tested by Thermo Fisher Scientific to ensure that the kits perform according to specifications and are free from contaminants that interfere with the assay performance and analysis.

The Quantifiler™ Duo DNA Quantification Kit can quantify DNA sample concentrations between 23 pg/L and 50 ng/L when using the reagents in the kit. Follow the protocols that are described in the *Quantifiler™ Duo DNA Quantification Kit User Guide* and run the samples on an Applied Biosystems™ 7500 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.

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. 4387913

Revision	Date	Description
F	22 October 2018	Updated branding and trademarks, no technical changes.
E	March 2012	Basis for this revision.

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. AmpliTaq Gold is a trademark of Roche Molecular Systems, Inc.