

Quantifiler® HP and Trio DNA Quantification Kits

Pub. No. 4485356 Rev. A

Note: For safety and biohazard guidelines, refer to the “Safety” section in the *Quantifiler® HP and Trio DNA Quantification Kits User Guide* (Pub. no. 4485354). For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

■ Product overview	1
■ Guidelines for calculating the standards dilution series	1
■ Prepare the DNA quantification standards	1
■ Prepare the reactions	2
■ Run the plate on the 7500 Real-Time PCR instrument	2
■ Analyze the experiment	3
■ View and export results	4
■ Limited Product Warranty	4

Product overview

This document describes the Quantifiler® HP DNA Quantification Kit (Cat. no. 4482911) and Quantifiler® Trio DNA Quantification Kit (Cat. no. 4482910). The Quantifiler® HP Kit is designed to quantify the total amount of amplifiable human DNA in a sample. The Quantifiler® Trio Kit is designed to simultaneously quantify the total amount of amplifiable human DNA and human male DNA in a sample.

Guidelines for calculating the standards dilution series

We recommend a ten-fold dilution series with 5 concentration points, for example, 0.005 ng/μL, 0.050 ng/μL, 0.500 ng/μL, 5.000 ng/μL, and 50.000 ng/μL. You can add an optional 100 ng/μL standard point if needed. However, you may see an increase in the IPC C_T for the 100 ng/μL standard. See the user guide for a standards dilution series example.

Use a minimum input volume of 10 μL DNA for dilutions (to ensure accuracy of manual pipetting).

Prepare the DNA quantification standards

When using Quantifiler® THP DNA Dilution Buffer, you can store the prepared DNA quantification standards in low-binding tubes for up to 2 weeks at 2 to 8°C. To prepare the DNA quantification standards dilution series:

1. Label five microcentrifuge tubes: Std. 1, Std. 2, Std. 3, and so on.
2. Dispense the required amount of diluent (Quantifiler® THP DNA Dilution Buffer) to each tube.
3. Prepare Std. 1:
 - a. Vortex the Quantifiler® THP DNA Standard 3 to 5 seconds.
 - b. Using a new pipette tip, add the appropriate volume of Quantifiler® THP DNA Standard for your dilution series to the tube for Std. 1.
 - c. Mix the dilution thoroughly.
4. Prepare Std. 2 through 5:
 - a. Using a new pipette tip, add the appropriate volume of the prepared standard to the tube for the next standard.
 - b. Mix the standard thoroughly.
 - c. Repeat steps a and b for each subsequent standard until you complete the dilution series.

Prepare the reactions

While preparing the reactions, keep the 96-well optical reaction plate or optical 8-tube strip in its base and do not place it directly on the bench top to protect it from scratches and particulate matter. To prepare the reactions:

1. Calculate the volume of each component needed to prepare the reactions, using the appropriate table below.

For the Quantifiler[®]HP and Trio DNA Quantification Kits:

Component	Volume Per Reaction (µL)
Primer Mix	8
Reaction Mix	10

Note: Include additional reactions in your calculations to provide excess volume for the loss that occurs during reagent transfers.

2. Prepare the reagents:
 - Thaw the Primer Mix completely, then vortex 3 to 5 seconds and centrifuge briefly before opening the tube.
 - Gently vortex the Reaction Mix before using.
3. Pipette the required volumes of components into an appropriately sized polypropylene tube.
4. Vortex the PCR mix 3 to 5 seconds, then centrifuge briefly.
5. Dispense 18 µL of the PCR mix into each reaction well or tube.
6. Add 2 µL of sample, standard, or control to the applicable wells or tubes.

Note: We recommend running duplicates of each sample of the DNA quantification standards for each reaction plate.
7. Seal the reaction plate with the Optical Adhesive Cover, or the strip tube with the optical 8-cap strip.
8. Centrifuge the plate at 3000 rpm for about 20 seconds in a tabletop centrifuge with plate holders to remove any bubbles.

Note: If a tabletop centrifuge with 96-well plate adapters is not available, visually inspect the plate for bubbles, and lightly tap the plate to remove bubbles in wells.

Run the plate on the 7500 Real-Time PCR instrument

Before you run the reactions, power on the computer, 7500 Real-Time PCR instrument, and software. Create an experiment for the run. To run the plate on the 7500 Real-Time PCR instrument:

1. Press the tray door to open it.
2. Load the plate into the plate holder in the instrument. Ensure that the plate is correctly aligned in the holder.
3. Load the 96-well optical plate with the notched A12 position at the top-right of the tray.
4. Close the tray door.
5. Apply pressure to the right side of the tray and at an angle to close the tray door.
6. In the HID Real-Time PCR Analysis Software, open the experiment that you set up for the run.
7. Click **Start Run**.

Analyze the experiment

Analyze a run after it is complete and reanalyze after you make any changes to the experiment, such as sample names.

To analyze an experiment:

1. Open the experiment for analysis.
2. Verify the analysis settings:
 - a. Click **Analysis Settings** in the upper-right corner of the window.
 - b. Click the **C_T Settings** tab.

The recommended C_T settings for each Quantifiler[®] kit are included in the experiment templates provided with the software. They are also listed in the *Quantifiler[®] HP and Trio DNA Quantification Kits User Guide*, and shown below. The recommended settings are those which were used in the validation experiments performed for each kit by Thermo Fisher Scientific.

- c. Verify that the settings are as shown below, then:
 - If the analysis settings are correct, click **Apply Analysis Settings**, or
 - If the analysis settings differ from those shown below, change them to match the settings, then click **Apply Analysis Settings**.

Large, small, and Y target threshold and baseline settings:

Target	Threshold	Baseline Start	Baseline End
T.IPC	0.1	3	15
T.Large Autosomal	0.2	3	15
T.Small Autosomal	0.2	3	15
T.Y	0.2	3	15

Cr Settings for T.Large Autosomal

Cr Settings to Use: Use Default Settings

Automatic Threshold

Threshold: 0.2

Automatic Baseline

Baseline Start Cycle: 3 End Cycle: 15

IPC target threshold and baseline settings:

Target	Threshold	Baseline Start	Baseline End
T.IPC	0.1	3	15
T.Large Autosomal	0.2	3	15
T.Small Autosomal	0.2	3	15
T.Y	0.2	3	15

Cr Settings for T.IPC

Cr Settings to Use: Use Default Settings

Automatic Threshold

Threshold: 0.1

Automatic Baseline

Baseline Start Cycle: 3 End Cycle: 15

For IPC threshold, 0.1
(see the note below)

Note: Quantifiler[®] HP and Trio Kits have been validated using the Manual Baseline method. Studies were also performed applying the Automatic Baseline method and the Manual Baseline method to evaluate potential differences between the methods for concentrations from 5 – 0.005 ng/μL. No statistically significant differences were observed within this range for C_T values generated using the Automatic Baseline and Manual analysis methods.

A value of 0.1 was used for the IPC Threshold during the developmental validation studies. Before using alternative baseline methods, (e.g. automatic) or thresholds, perform the appropriate internal validation studies.

3. Click **Analyze**.

View and export results

For information about how to view and export your results, see the *Quantifiler[®] HP and Trio DNA Quantification Kits User Guide* (Pub. no. 4485354).

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.

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

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

DISCLAIMER: 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.

© 2014 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com

29 April 2014

