



Brilliant III Ultra-Fast QRT-PCR Master Mix

Quick Reference Guide for the LightCycler® 480 Real-Time PCR System

This quick reference guide provides an optimized protocol for using Agilent's Brilliant III Ultra-Fast QRT-PCR Master Mix with the LightCycler 480 Real-Time PCR System from Roche. For detailed instructions, refer to the full product manual.

Prepare the Reactions

- 1 Prepare the experimental reactions by combining the components of the reagent mixture in the order listed in the table below. Prepare a single reagent mixture for replicate reactions (plus at least one reaction volume excess) using multiples of each component. *Keep reagent mixture on ice.*

| Reagent Mixture |
|---|
| Nuclease-free PCR-grade water to bring final volume to 20 μ l (including RNA) |
| 10 μ l of 2 \times QRT-PCR Master Mix |
| x μ l of experimental probe at optimized concentration (100–600 nM) |
| x μ l of upstream primer at optimized concentration (200–600 nM) |
| x μ l of downstream primer at optimized concentration (200–600 nM) |
| 0.2 μ l of 100 mM DTT |
| 1 μ l of RT/RNase Block |

- 2 Gently mix the reagent mixture without creating bubbles, then distribute the mixture to the experimental reaction tubes. *Keep the reactions on ice.*
- 3 Add x μ l of experimental RNA to each reaction to bring the final reaction volume to 20 μ l. The table below lists a suggested quantity range for different RNA templates.

| RNA | Quantity per reaction |
|-----------|-----------------------|
| Total RNA | 0.1 pg – 100 ng |
| mRNA | 0.1 pg – 1 ng |

- 4 Mix the reactions without creating bubbles, then centrifuge briefly.



Set Up the QPCR Plate and Thermal Profile

- 1 From the main window in the LightCycler 480 software, click **Sample Editor** on the module bar to open the *Sample Editor* module. Enter sample information for your experiment as needed.
- 2 Click **Experiment** on the module bar to open the *Run* module.
- 3 From the **Run Protocol** tab, enter a reaction volume of 20 μ l.
- 4 Set the **Detection Format** to *Mono Color Hydrolysis Probe* or *Multi Color Hydrolysis Probe* as appropriate for your experiment.
- 5 Set up the PCR program to run the cycling protocol below:

| Program Name | Cycles | Analysis Mode | Acquisition Mode | Ramp Rate ($^{\circ}$ C/s) | Hold Time | Temperature |
|-----------------------|--------|----------------|------------------|-----------------------------|------------|-----------------|
| Reverse transcription | 1 | None | None | 4.4 | 10 minutes | 50 $^{\circ}$ C |
| Denaturation | 1 | None | None | 4.4 | 3 minutes | 95 $^{\circ}$ C |
| Amplification | 45 | Quantification | None | 4.4 | 5 seconds | 95 $^{\circ}$ C |
| | | | Single | 2.2 | 10 seconds | 60 $^{\circ}$ C |
| Cooling | 1 | None | None | 2.2 | 30 seconds | 40 $^{\circ}$ C |

Run the PCR Program

- 1 Place the reactions in the LightCycler 480 instrument.
- 2 From the **Run Protocol** or **Data** tab, click **Start Run**.

Analyze Data

- 1 Analyze the results of the run as needed for your experiment.

Notice to Purchaser

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

LightCycler[®] is a registered trademark of Roche.

Product Information

Catalog #600884, 400 reactions
Catalog #600885, 4000 reactions

Ordering Information

By phone (US and Canada*): 800-227-9770
On the web: www.agilent.com/genomics

Technical Services

By phone (US and Canada*): 800-227-9770
By email: techservices@agilent.com

*For other countries, please contact your local sales representative at www.agilent.com/genomics/contactus