



# Brilliant III Ultra-Fast SYBR<sup>®</sup> Green QPCR Master Mix with High ROX

## Quick Reference Guide for the ABI 7900HT Fast Real-Time PCR System

*This quick reference guide provides an optimized protocol for using Agilent's Brilliant III Ultra-Fast SYBR<sup>®</sup> Green QPCR Master Mix with High ROX (Catalog #600889) with the 7900HT Fast Real-Time PCR System from Life Technologies. 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.

Reagent Mixture
Nuclease-free PCR-grade water to bring final volume to 20 $\mu$ L (including DNA)
10 $\mu$ L of 2 $\times$ SYBR Green QPCR Master Mix
$x$ $\mu$ L of upstream primer at optimized concentration (200–500 nM)
$x$ $\mu$ L of downstream primer at optimized concentration (200–500 nM)

- 2 Gently mix the reagent mixture without creating bubbles, then distribute the mixture to the experimental reaction tubes.
- 3 Add  $x$   $\mu$ L of experimental DNA to each reaction to bring the final reaction volume to 20  $\mu$ L. The table below lists a suggested quantity range for different DNA templates.

DNA	Quantity per reaction
Genomic DNA	5 $\mu$ g – 50 ng
cDNA	0.5 $\mu$ g – 100 ng*

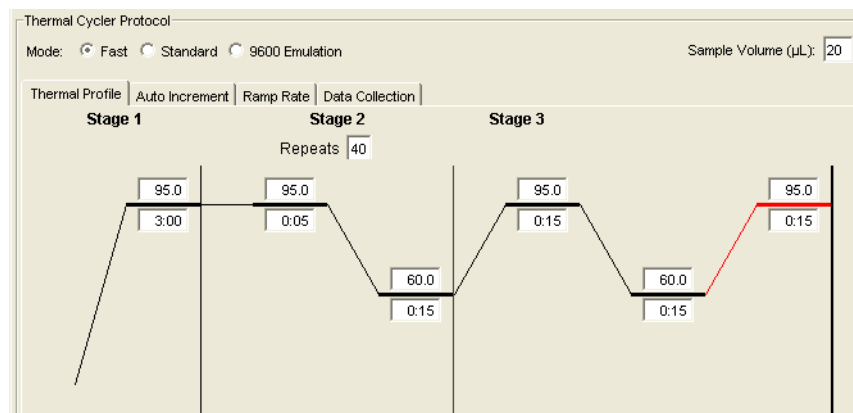
\*Refers to RNA input amount during cDNA synthesis

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



## Set Up the QPCR Plate and Thermal Profile

- 1 From the SDS software, click **File > New** to open the Plate Document Wizard.
- 2 Enter the appropriate assay and well information for a new experiment.
- 3 Click **OK**. The Wizard will close and the plate document will appear in the main software window.
- 4 Click **Add Detector**, and select the correct reporter for the assay. Click **Copy to Plate Document**, then click **Done**.
- 5 Highlight the wells that will contain samples and check the selected reporter dye.
- 6 On the Instrument/Thermal Profile tab, enter a sample volume of 20  $\mu\text{L}$  and select the *Fast* run mode. Adjust the thermal cycling conditions according to the image below, and set the instrument to report fluorescence during the 60°C step of each cycle.



## Run the PCR Program

- 1 Place the reactions in the 7900HT instrument.
- 2 On the Instrument/Real Time tab, click **Start Run**.

## Analyze Data

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

**Endnote:** SYBR<sup>®</sup> Green is a registered trademark of Molecular Probes, Inc.

### Product Information

Catalog #600889, 400 reactions

### Ordering Information

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

### Technical Services

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

\*For other countries, please contact your local sales representative at [www.agilent.com/genomics/contactus](http://www.agilent.com/genomics/contactus)

Manual Part Number 5990-7251, Revision B0  
For Research Use Only. Not for use in diagnostic procedures.

©Agilent Technologies, Inc. 2013, 2016