



Brilliant III Ultra-Fast SYBR® Green QPCR 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 SYBR® Green QPCR 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.

Reagent Mixture
Nuclease-free PCR-grade water to bring final volume to 20 µl (including DNA)
10 µl of 2× SYBR Green QPCR Master Mix
x µl of upstream primer at optimized concentration (200–500 nM)
x µl 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 µl of experimental DNA to each reaction to bring the final reaction volume to 20 µl. The table below lists a suggested quantity range for different DNA templates.

DNA	Quantity per reaction
Genomic DNA	5 pg – 50 ng
cDNA	0.5 pg – 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 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 µl.
- 4 Set the **Detection Format** to *SYBR Green I/HRM Dye*.
- 5 Set up the PCR program to run the cycling protocol below:

Program Name	Cycles	Analysis Mode	Acquisition Mode	Ramp Rate (°C/s)	Hold Time	Temperature
Pre-incubation	1	None	None	4.4	3 minutes	95°C
Amplification	45	Quantification	None	4.4	5 seconds	95°C
			Single	2.2	10 seconds	60°C
Melting curve	1	Melting curves	None	4.4	5 seconds	95°C
			None	2.2	1 minute	65°C
			Continuous (5 acquisitions/second)	0.11	—	97°C
Cooling	1	None	None	2.2	30 seconds	40°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.

Notices to Purchaser

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. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

SYBR® is licensed for research and development only under patents and patent applications owned by Invitrogen Corporation.

LightCycler® is a registered trademark of Roche.

SYBR® is a registered trademark of Molecular Probes, Inc.

Product Information

Catalog #600882, 400 reactions
Catalog #600883, 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

Manual Part Number 5990-7210, Revision B0
For Research Use Only. Not for use in diagnostic procedures.

©Agilent Technologies, Inc. 2015
5990-6144EN