



# Brilliant III Ultra-Fast SYBR® Green QPCR Master Mix with Low ROX

## Quick Reference Guide for the Agilent Mx3000P/Mx3005P QPCR Systems

*This quick reference guide provides an optimized protocol for using Agilent's Brilliant III Ultra-Fast SYBR® Green QPCR Master Mix with Low ROX (Catalog #600892) with the Mx3000P and Mx3005P QPCR Systems. 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*

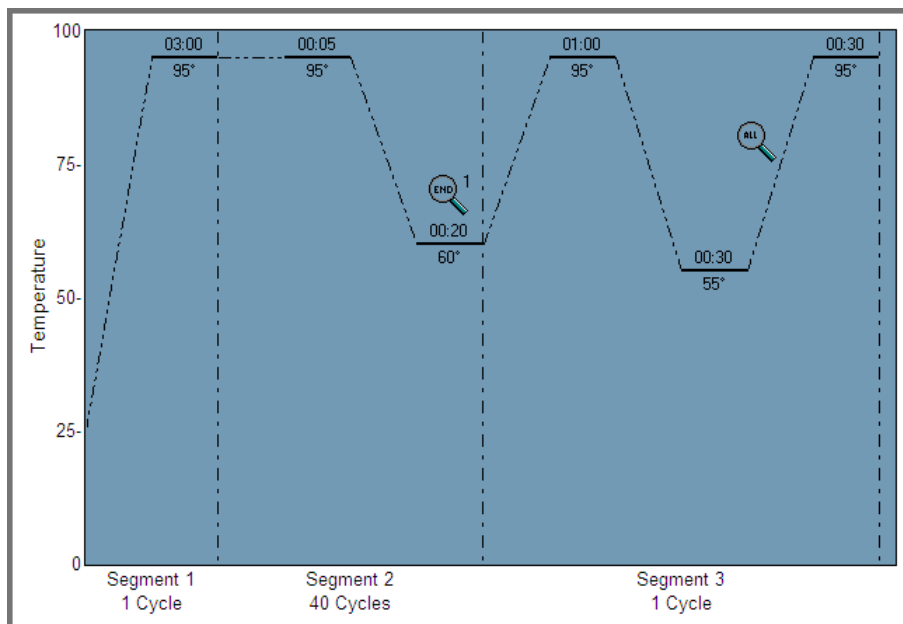
\*Refers to RNA input amount during cDNA synthesis

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



## Set Up the QPCR Plate and Thermal Profile

- 1 Complete the **Plate Setup** screen for a new experiment as needed, including assigning well types and assay information.
- 2 On the **Thermal Profile Setup** screen, set the **Thermal Profile Design** selection to **Standard**.
  - Under **Pre-Melt/RT Segment**, click **1 Plateau**.
  - Under **Amplification Segment**, click **Fast 2 Step**.
  - Under **Dissociation/Melt Segment**, click **Dissociation/Melt**.
- 3 Adjust the thermal profile according to the image below. The profile includes a 5-second denaturation step. Note that some assays may require a denaturation of up to 20 seconds. The exact denaturation time needs to be optimized for each target.



## Run the PCR Program

- 1 Place the reactions in the Mx3000P/Mx3005P instrument.
- 2 On the **Run** screen, click **Start Run**.

## Analyze Data

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

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

### Product Information

Catalog #600892, 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)