



Brilliant III Ultra-Fast SYBR[®] Green QRT-PCR Master Mix

Quick Reference Guide for the Stratagene Mx3000P/Mx3005P QPCR Systems

This quick reference guide provides an optimized protocol for using the Stratagene Brilliant III Ultra-Fast SYBR[®] Green QRT-PCR Master Mix with the Stratagene Mx3000P and Mx3005P QPCR Systems from Agilent. For detailed instructions, refer to the full product manual.

Prepare the Reactions

- 1 Dilute the reference dye 1:500 using nuclease-free PCR-grade water.
- 2 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 the 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 SYBR Green QRT-PCR Master Mix
x μ l of upstream primer at optimized concentration (150–500 nM)
x μ l of downstream primer at optimized concentration (150–500 nM)
0.3 μ l of diluted reference dye
0.2 μ l of 100 mM DTT
1 μ l of RT/RNase Block

- 3 Gently mix the reagent mixture without creating bubbles, then distribute the mixture to the experimental reaction tubes. *Keep the reactions on ice.*
- 4 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.

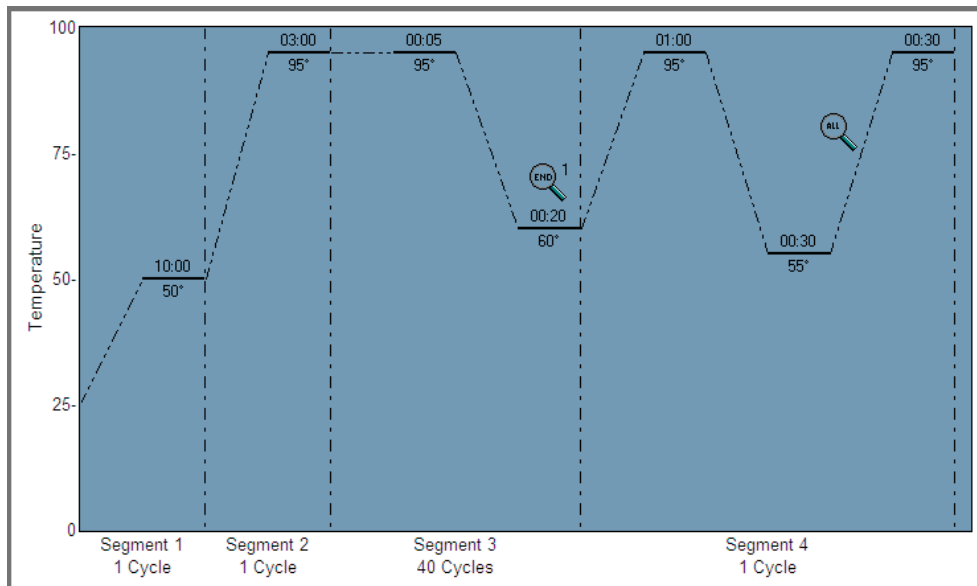
RNA	Quantity per reaction
Total RNA	0.1 pg – 100 ng
mRNA	0.1 pg – 1 ng

- 5 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 **2 Plateaus**.
 - 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.

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.

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

Product Information

Catalog #600886, 400 reactions
Catalog #600887, 4000 reactions

Ordering Information

By phone (US only*): 800-424-5444, x3
On the web: www.genomics.agilent.com

Technical Services

By phone (US only*): 800-894-1304, x2
By email: techservices@agilent.com

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

Manual Part Number 5990-7207, Revision A

For Research Use Only. Not for use in diagnostic procedures.

©Agilent Technologies, Inc. 2010

Printed June 1, 2010

5990-5997EN