



# Brilliant III Ultra-Fast SYBR<sup>®</sup> Green QRT-PCR Master Mix

## Quick Reference Guide for the ABI 7500 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 QRT-PCR Master Mix with the 7500 Fast Real-Time PCR System from Applied Biosystems. 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 $\mu$ l (including RNA)
10 $\mu$ l of 2 $\times$ SYBR Green QRT-PCR Master Mix
x $\mu$ l of upstream primer at optimized concentration (150–500 nM)
x $\mu$ l of downstream primer at optimized concentration (150–500 nM)
0.3 $\mu$ l of diluted reference dye
0.2 $\mu$ l of 100 mM DTT
1 $\mu$ 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  $\mu$ l of experimental RNA to each reaction to bring the final reaction volume to 20  $\mu$ 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.



