

# TAQuest™ qPCR Master Mix for TaqMan Probes \*No ROX\*

Catalog number: 17282, 17283 Unit size: 1 mL, 5 mL

Component	Storage	Amount (Cat No. 17282)	Amount (Cat No. 17283)
TAQuest™ qPCR Master Mix for TaqMan Probes	Freeze (< -15 °C), Minimize light exposure	1 mL	5 mL
*No ROX*			

#### **OVERVIEW**

TAQuest™ qPCR Master Mix for TaqMan Probes is a ready-to-use 2X solution optimized for qPCR and 2-step RT-qPCR compatible with TaqMan gene expression assays. The master mix provides all of the essential components including our proprietary TAQuest™ Hot Start Taq DNA Polymerase enzyme and dNTPs in an optimized PCR buffer, except the template, primers and probes. The Hot Start Taq DNA polymerase allows you to set up a PCR reaction at room temperature, thus minimizing non-specific product formation. The optimized composition ensures PCR specificity and sensitivity with all sample types such as genomic, plasmid, viral and cDNA templates. TAQuest™ qPCR Master Mix for TaqMan Probes has been designed to be used for duplex reactions using internal positive controls. This master mix does not contain a ROX reference dye.

#### **KEY PARAMETERS**

#### qPCR

Instrument specification(s)

Filter based on probes

## SAMPLE EXPERIMENTAL PROTOCOL

The following protocol can be used as a guideline.

**Note** Thaw the TAQuest™ qPCR Master Mix for TaqMan Probes \*No ROX\* at room temperature. Vortex qPCR Master Mix thoroughly before use.

- 1. Prepare one of the following reaction mixes as indicated in Table 1.
- Carefully mix the reagents with a gentle vortex followed by a brief centrifuge.
- Set up the plate in the qPCR instrument and run as indicated in Table
  2.

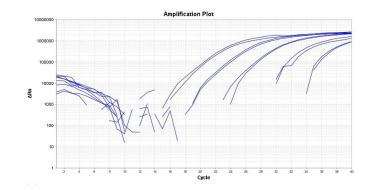
Table 1. Reagents composition per well for each reaction

Components	Volume (25 µ L/reaction)	Volume (50 μ L/reaction)	Final Conc.
TAQuest™ qPCR Master Mix for TaqMan Probes *No ROX*	12.5 μL	25 μL	1X
Upstream primer, 10 µM	0.25-2.5 μL	0.5-5.0 μL	0.1-1.0 μM
Downstream primer, 10 μΜ	0.25-2.5 μL	0.5-5.0 μL	0.1-1.0 μM
TaqMan Probes, 10 μM	0.25-0.625 µL	0.5-1.25 μL	100-250 nM
DNA template	1-5 µL	1-5 µL	Optimized conc.
Nuclease-Free Water to	25 μL	50 μL	

Table 2. Thermal cycling parameters

Parameter	Polymerase Activation	PCR (30-40 cycles)		
	Hold	Denature	Anneal	Extend
Temperature	95 °C	95 °C	55-65 °C	68-72 °C
Time (m:ss)	0:20	0:30	1:00	1:00

# EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Amplification plot for a dilution series of HeLa cells cDNA amplified in replicate reactions to detect GAPDH using TAQuest™ qPCR Master Mix for TaqMan Probes \*No ROX\*.

## **DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.