

# TaqMan® Small RNA Assays

## TaqMan® MicroRNA Assays, TaqMan® siRNA Assays, and Custom TaqMan® Small RNA Assays

**Note:** For safety and biohazard guidelines, refer to the “Safety” section in the *TaqMan® Small RNA Assays Protocol* (Part no. 4364031). For all chemicals, read the SDS and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

### Perform reverse transcription (RT)

- 1** Prepare the total RNA
  - a. Isolate the total RNA.  
Use a method that preserves small RNAs. To prevent the loss of the longer control transcripts (snoRNAs), we recommend that you do *not* perform size fractionation.
  - b. Quantify the amount of total RNA in your sample.

- 2** Prepare the RT master mix
  - a. Allow the TaqMan® MicroRNA Reverse Transcription Kit components to thaw on ice.
  - b. In a polypropylene tube, prepare the RT master mix on ice by scaling the volumes listed below to the desired number of RT reactions. We recommend adding 10–20% overage to account for pipetting losses.

Component	Master mix volume per 15-µL reaction†
100mM dNTPs (with dTTP)	0.15 µL
MultiScribe™ Reverse Transcriptase, 50 U/µL	1.00 µL
10X Reverse Transcription Buffer	1.50 µL
RNase Inhibitor, 20 U/µL	0.19 µL
Nuclease-free water	4.16 µL
<b>Total volume</b>	<b>7.00 µL</b>

† Each 15-µL RT reaction consists of 7 µL master mix, 3 µL of 5X RT primer, and 5 µL RNA sample.

- 3** Prepare the RT reaction
  - a. Thaw the 5X RT primer and RNA template on ice. Before use, vortex the RT primer tubes to mix, then centrifuge briefly.
  - b. If you are performing quantitation of:
    - Ambion *Silencer*® Select siRNAs, go to [step c](#).
    - All other templates, go to [step d](#).
  - c. Mix gently. Centrifuge to bring solution to the bottom of the tube.
  - d. Place the RT master mix on ice until you prepare the RNA reaction.

### 3 Prepare the RT reaction (continued)

- c. If you are performing Ambion *Silencer*® Select siRNA quantitation, denature and prepare the double-stranded template:
1. For each 15- $\mu$ L RT reaction, combine 3  $\mu$ L of 5X RT primer and 5  $\mu$ L of double-stranded template in a 0.2-mL reaction tube or a well of a 96-well reaction plate.
  2. Incubate the tube or plate at 85°C for 5 minutes, then at 60°C for 5 minutes.
  3. Place the denatured template on ice.
  4. For each 15- $\mu$ L RT reaction, combine RT master mix with denatured RNA and RT primer in the ratio of:  
7  $\mu$ L RT master mix : 8  $\mu$ L denatured RNA and RT primer (1–10 ng of RNA per reaction)
  5. Go to [step e](#).
- d. If you are preparing single-stranded RNA, prepare the total RNA template:
1. For each 15  $\mu$ L RT reaction, combine RT master mix with 1–10 ng total RNA in the ratio of:  
7  $\mu$ L RT master mix : 5  $\mu$ L total RNA
  2. Mix gently, then centrifuge briefly to bring the solution to the bottom of the tube.
  3. Transfer 12.0  $\mu$ L of the RT master mix-total RNA mixture to a 0.2-mL reaction tube or a well of a 96-well reaction plate.
  4. Add 3  $\mu$ L of 5- $\mu$ L RT primer from each assay set to the corresponding RT reaction tube or well.
  5. Go to [step e](#).
- e. Seal the reaction tube or reaction plate, then mix gently. Centrifuge briefly.
- f. Incubate the reactions on ice for 5 minutes, then keep on ice until you load the thermal cycler.

### 4 Perform reverse transcription

Load the reaction tubes or plate into a thermal cycler, then perform a run using the following conditions:

- Mode: Standard
- Reaction volume: 15  $\mu$ L
- Thermal cycling conditions:

Step	Time	Temperature
Hold	30 minutes	16°C
Hold	30 minutes	42°C
Hold	5 minutes	85°C
Hold	$\infty$	4°C

## Quantitative PCR (qPCR) amplification

### 1 Prepare the qPCR reactions

- Place the components on ice. Gently invert the tubes to mix, then return them to the ice.
- Calculate the volumes needed based on the number of reactions and a reaction volume of 20  $\mu\text{L}$ .

**Note:** We recommend performing three replicates of each reaction. Include excess volume in your calculations to account for loss that occurs during reagent transfers.

Component	Volume ( $\mu\text{L}$ )	
	20- $\mu\text{L}$ reaction	3 replicates + 20% excess
TaqMan® Universal PCR Master Mix II, no UNG <sup>†</sup>	10.00 $\mu\text{L}$	36.00 $\mu\text{L}$
Nuclease-free water	7.67 $\mu\text{L}$	27.61 $\mu\text{L}$
TaqMan® Small RNA Assay (20X)	1.00 $\mu\text{L}$	3.60 $\mu\text{L}$
Product from RT reaction	1.33 <sup>‡</sup> $\mu\text{L}$	4.80 $\mu\text{L}$
<b>Total volume</b>	<b>20.00 <math>\mu\text{L}</math></b>	<b>72.01 <math>\mu\text{L}</math></b>

<sup>†</sup> TaqMan Universal PCR Master Mix II with UNG is compatible with TaqMan® Small RNA Assays.

<sup>‡</sup> Maximum volume of RT product that can be added to each reaction.

- Combine reaction components in a microcentrifuge tube.
- Mix gently by inversion, then centrifuge the tube or reaction plate briefly.

### 2 Prepare the qPCR reaction plate

- Transfer 20  $\mu\text{L}$  of the PCR reaction into wells of a reaction plate.
- Seal the plate with optical adhesive film or optical caps, then briefly centrifuge the reaction plate.

### 3 Run the PCR reaction plate

- Create an experiment or plate document using the following parameters:
  - Mode: Standard
  - Reaction volume: 20  $\mu\text{L}$
  - Thermal cycling conditions:

Step	Optional AmpErase® UNG activity <sup>†</sup>	Enzyme Activation	PCR	
	HOLD	HOLD	CYCLE (40 cycles)	
			Denature	Anneal/extend
Temperature	50°C	95°C	95°C	60°C
Time	2 minutes	10 minutes	15 seconds	60 seconds

<sup>†</sup> Not needed when UNG is not in the reaction.

- Load the reaction plate into the real-time PCR instrument.
- Start the run.

### 4 Analyze the experiment

Refer to the getting started guides for your real-time PCR system to analyze the experiment. The general process for analyzing the data from gene expression assays involves the following procedures:

- View the amplification plots.
- Set the baseline and threshold values.

---

**For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.**

NOTICE TO PURCHASER: PLEASE REFER TO THE TAQMAN® SMALL RNA ASSAYS SUBMISSION GUIDELINES PROTOCOL FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION

Trademarks of Life Technologies Corporation and its affiliated companies: AB (Design)®, Applied Biosystems®, MultiScribe™, Silencer®, TaqMan, AmpErase and AmpliTaq Gold are registered trademarks of Roche Molecular Systems, Inc. TRI Reagent is a registered trademark of Molecular Research Center, Inc.

All other trademarks are the sole property of their respective owners.

© 2011 Life Technologies Corporation Corporation. All rights reserved.

Part Number 4412551 Rev. C 01/2011



**Headquarters**  
5791 Van Allen Way | Carlsbad, CA 92008 USA  
Phone 760.603.7200  
[www.lifetechnologies.com](http://www.lifetechnologies.com)

**Technical Resources and Support**  
For the latest technical resources and support information  
for all locations, please refer to our Web site at  
[www.appliedbiosystems.com](http://www.appliedbiosystems.com)