

I. Introduction

Lyo-Ready One Step PrimeScript III RT-qPCR Mix (Cat. No. 638345) enables the user to perform accurate probe-based RT-qPCR on purified RNA samples using a lyophilized master mix. Primers and probes can be added to the mix prior to lyophilization. Once dried, the master mix allows for flexibility in handling larger sample volumes, in addition to the convenience of room-temperature storage and transport.

II. Required Materials

This protocol applies to the following Takara Bio products:

- Lyo-Ready One Step PrimeScript III RT-qPCR Mix*

*Store at –20°C.

Additional materials required:

- Lyophilizer (e.g., SP VirTis Genesis Pilot Freeze Dryer, SP VirTis Ultra Pilot-Production Freeze Dryer (ATS Automation Tooling Systems, Inc.), etc.)
- Primers and probes
- Micropipette and tips (with hydrophobic filters)
- Vortex mixer
- Benchtop centrifuge for tubes or plates
- 1.5 ml Eppendorf tubes, 200 µl PCR tubes, or 200 µl PCR plates for sample preparation
- Tubes or plates for real-time PCR with optical seals or caps
- A real-time PCR machine (e.g., CFX96 Real-Time PCR Detection System (Bio-Rad), QuantStudio 3 or 5 (Thermo Fisher Scientific))

III. Protocol

A. Lyophilization

The master mix can be lyophilized in concentrations from 1X to 2X and can include primers and probes (i.e., assay). Lyophilization recipes will vary based on individual needs. Optimization will be required.

- Total cycle time and drying times will vary; 15–24+ hours can be expected.
- Long-term storage at ambient temperatures requires that the lyophilized product be packaged in a heat-sealed pouch with silica desiccant while at low relative humidity conditions.

Example:

1. Dispense 10 µl Lyo-Ready One Step PrimeScript III RT-qPCR Mix into each tube of an 8-well strip in a 96-well rack.
2. With the lids open on the tubes, place racks into a lyophilizer.
3. Freeze at –70°C for 4 hr.
4. Dry at –30°C, 150 mTorr until primary drying is complete (≥6.5 hr).
5. Perform secondary drying at 20–25°C.

B. RT-qPCR Reaction Mix Preparation

1. Prepare the reaction mix for the RT-qPCR reaction as illustrated below (optimization is recommended). After preparation, vortex the mixture, centrifuge, and immediately place the tube/plate into the thermal cycler.

NOTE: Adding up to 10% DMSO can improve performance with some assays (optimization required).

Examples:

- **RT-qPCR reaction mixture (1 rxn) for cakes lyophilized with an assay**

1 cake	Lyophilized One Step PrimeScript III RT-qPCR Mix+Primers/Probes
18 µl	RNase Free H ₂ O
2 µl	RNA sample*
20 µl	Total volume

- **RT-qPCR reaction mixture (1 rxn) for cakes without an assay**

1 cake	Lyophilized One Step PrimeScript III RT-qPCR Mix
1 µl	20X primer/probe mix (provided by user)
17 µl	RNase Free H ₂ O
2 µl	RNA sample*
20 µl	Total volume

*The RNA range will be dependent on your individual assay and testing conditions. Please experiment to find the volume optimal for your experimental parameters.

2. Vortex until cake is completely solubilized then spin down.
3. Wait 5 min, vortex, and spin down again.

C. RT-qPCR Reaction Protocol

1. Run the standard cycling protocol according to the instruction manual of your real-time qPCR instrument.

NOTE: The PCR program below was used for the CDC’s SARS-CoV-2 multiplex assay of three targets (SARS-CoV-2 N1 and N2 targets and Human RNaseP target) and high-quality results were obtained. Further optimization may be required for your experiment.

52°C	5 min	Reverse transcription
95°C	2 min	Initial denaturation
40 cycles:		
95°C	10 sec	Denaturation Anneal, extend, and capture
60°C	30 sec	

2. After the reaction is complete, check the amplification curve. Confirm that the analytical parameters are appropriate and that the Cq value has been calculated.

NOTE: If the default setting analysis does not work, perform manual analysis per the instrument’s instruction manual.

Contact Us	
Customer Service/Ordering	Technical Support
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)
web: takarabio.com/service	web: takarabio.com/support
e-mail: ordersUS@takarabio.com	e-mail: technical_support@takarabio.com

Notice to Purchaser

Our products are to be used for **Research Use Only**. They may not be used for any other purpose, including, but not limited to, use in humans, therapeutic or diagnostic use, or commercial use of any kind. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without our prior written approval.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

© 2023 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

This document has been reviewed and approved by the Quality Department.