

## I. Introduction

**Lyo-Ready PrimePath™ Probe qPCR Mix, GPR** (Cat. No. 638352) enables the user to perform accurate probe-based qPCR on purified DNA 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 PrimePath Probe qPCR Mix, GPR (2X)\*

\*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 and at low relative humidity conditions.

### Example:

1. Dispense 10 µl Lyo-Ready PrimePath Probe qPCR Mix, GPR 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 ( $\geq 6.5$  hr).
5. Perform secondary drying at 20–25°C.

**SAFE STOPPING POINT:** Lyophilized cakes can be stored long-term following the parameters above.

## B. qPCR Reaction Mix Preparation

1. Prepare the reaction mix for the qPCR reaction as illustrated below (optimization is recommended).

### Examples:

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

1 cake	Lyophilized PrimePath Probe qPCR Mix + primers/probes
18 µl	DNase/RNase Free H <sub>2</sub> O
2 µl	DNA sample*
<b>20 µl</b>	<b>Total volume</b>

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

1 cake	Lyophilized PrimePath Probe qPCR Mix
1 µl	20X primer/probe mix (provided by user)
17 µl	DNase/RNase Free H <sub>2</sub> O
2 µl	DNA sample*
<b>20 µl</b>	<b>Total volume</b>

\*The DNA range will be dependent on your individual assay and testing conditions. You will need to optimize the volume for your experimental parameters.

2. Vortex until cake is completely dissolved then spin down.
3. Wait 5 min, then repeat Step 2.

**NOT A SAFE STOPPING POINT:** Proceed to the qPCR reaction protocol (Section C).

## C. qPCR Reaction Protocol

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

The example PCR program below was used for detecting a multiplex assay of five targets (*Chlamydia trachomatis*, *Neisseria gonorrhoeae*, *Trichomonas vaginalis*, an spike-in control, and Human RNaseP target) and high-quality results were obtained.

**NOTE:** Further optimization may be required for your experiment.

### Example:

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 C<sub>q</sub> 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: <a href="http://takarabio.com/service">takarabio.com/service</a>	web: <a href="http://takarabio.com/support">takarabio.com/support</a>
e-mail: <a href="mailto:ordersUS@takarabio.com">ordersUS@takarabio.com</a>	e-mail: <a href="mailto:technical_support@takarabio.com">technical_support@takarabio.com</a>

## Notice to Purchaser

This product is a general-purpose reagent intended **For Laboratory Use**. Outside of the United States, this product is intended for research use only unless otherwise stated. This product is not intended for a specific application or made for any particular clinical use. The performance characteristics of this product have not been fully established. It is the user's responsibility to validate the performance of the product, and any component thereof, for any particular use. Resale or transfer of this product, any component thereof, or any substance produced through use of this product, or any component thereof, to any third party is expressly forbidden. To obtain additional rights, please contact your local office or go to our website for additional information.

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](http://takarabio.com). It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

© 2024 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](http://takarabio.com).

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