

Power SYBR[®] Green PCR Master Mix: Executed on Bio-Rad iQ5 Real-Time PCR Detection System

For safety and biohazard guidelines, refer to the “Safety” section in the *Power SYBR[®] Green PCR Master Mix and RT-PCR Protocol* (PN 4367218). Read the MSDS and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

This quick reference card provides simplified procedures for using the *Power SYBR Green PCR Master Mix* for real-time PCR assays on the iQ5 Real-Time PCR Detection System. The *Power SYBR Green PCR Master Mix and RT-PCR Protocol* (PN 4367218) provides detailed real-time PCR and RT-PCR procedures and ordering information for the *Power SYBR Green PCR Master Mix*.

1

Prepare the PCR reagents mix

- Allow the *Power SYBR Green PCR Master Mix* to thaw completely. Mix gently.
- In a polypropylene tube, prepare the PCR reagents mix by scaling the volumes listed below to the desired number of PCR reactions.

Note: Include extra volume to account for pipetting losses.

Reaction Component	20 μ L/Reaction	Final Concentration
<i>Power SYBR Green PCR Master Mix (2X)</i>	10	1X
Reverse Primer	Variable	50 to 300 nM
Forward Primer	Variable	50 to 300 nM
Template	Variable	1 to 100 ng
Nuclease-free water	Variable	-

- Mix gently. Do not vortex. Centrifuge briefly and then prepare the PCR reaction plate.

2

Set up the plate document

See your instrument user’s manual for detailed instructions on how to configure the plate documents.

The thermal-cycling condition for the *Power SYBR Green PCR Master Mix* are described in the table below:

Step	AmpliTaq Gold [®] Enzyme Activation (*)	PCR	
	Hold	Cycles (40 cycles)	
		Denature	Anneal/Extend (**)
Time	10 min	15 sec	60 sec
Temp (°C)	95	95	60

Another option is to download the template “ABPowerSYBRThermalProtocolIQ5” from www.appliedbiosystems.com. To use this template, open the software and follow the steps below:

- In the “set-up” tab, “protocol”, choose the protocol file named “ABPowerSYBRThermalProtocolIQ5”.
- In the “plate” tab, click “edit” button to select the dye*** and set-up the plate. Click “Save and Exit plate editing”.

Note:

* Never decrease the temperature or time for the AmpliTaq Gold[®] Enzyme Activation.

** Before setup, check your primer annealing temperature. If primer T_m is < 60°C, we recommend using a 3-step protocol.

*** ROX[™] is included in the *Power SYBR Green PCR Master Mix*, but will not be used for fluorescence normalization in the Bio-Rad iQ5 Manager software.

3	Run the PCR reaction plate	Load the reaction plate into the instrument, then click “Start Run” tab to start the run. See your instrument user’s manual for detailed instructions on how to load and run the plate.
4	Analyze the results	Data Analysis varies depending on the instrument. See the <i>Power SYBR</i> [®] Green PCR Master Mix and RT-PCR Protocol (PN 4367218) and your instrument user’s manual for detailed instructions on how to analyze the data.

Power SYBR[®] Green PCR Master Mix Products

Item	Part Number	Contents
<i>Power SYBR</i> Green PCR Master Mix ± <ul style="list-style-type: none"> • Mini-Pack • 1-Pack • 2-Pack • 5-Pack • 10-Pack • Bulk Pack 	<ul style="list-style-type: none"> • 4368577 • 4367659 • 4368706 • 4368702 • 4368708 • 4367660 	<ul style="list-style-type: none"> • 1 x 1 mL tube (40 reactions) • 1 x 5 mL tube (200 reactions) • 2 x 5 mL tube (400 reactions) • 5 x 5 mL tube (1000 reactions) • 10 x 5 mL tube (2000 reactions) • 1 x 50 mL tube (2000 reactions)
<ul style="list-style-type: none"> • <i>Power SYBR</i> Green RT-PCR Reagents Kit 	<ul style="list-style-type: none"> • 4368711 	<ul style="list-style-type: none"> • <i>Power SYBR</i> Green RT-PCR Reagents Kit (200 x 50 µL reactions) • TaqMan[®] Reverse Transcription Reagents (200 x 10 µL reactions)
Related Documentation <ul style="list-style-type: none"> • Protocol • Quick Reference Card 	<ul style="list-style-type: none"> • 4367218 • 4367219 	<ul style="list-style-type: none"> • 1 protocol • 1 card

± Based on 50 µL reaction volume.

© 2008 Applied Biosystems Inc. All rights reserved. **For Research Use Only. Not for use in diagnostic procedures.**
 NOTICE TO PURCHASER: Please refer to the *Power SYBR* Green PCR Master Mix and RT-PCR Protocol (PN 4367218) user’s manual FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION.
 Applied Biosystems and AB (Design) are registered trademarks and ROX is a trademark of Applied Biosystems Inc. or its subsidiaries in the U.S. and/or certain other countries. iQ5 is a trademark of Bio-Rad, Inc. SYBR is a registered trademark of Molecular Probes, Inc. AmpliTaq Gold and TaqMan are registered trademarks of Roche Molecular Systems, Inc. All other trademarks are the sole property of their respective owners.