



## Do you have all the pieces of the puzzle?

It's never been easier to get everything you need for your real-time PCR experiments.

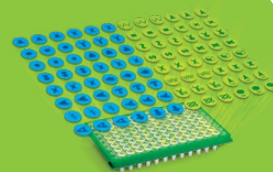
### 1 Cell Lysis and RNA Isolation

Aurum Total RNA Kits produce high-quality, DNA-free total RNA from various starting materials.



### 2 Target Selection

PrimePCR Assays and Panels are expertly designed PCR primer and probe assays that have been experimentally validated for guaranteed performance.



### 3 Reaction Setup

**Supermixes** — patented\* Sso7d fusion polymerase technology provides robust performance.

**Plastics** — patented Hard-Shell PCR Plates provide superior stability and flatness.



### 4 Detection

CFX Real-Time PCR Systems are designed to bring outstanding thermal performance and sensitive optical detection to all of your real-time PCR applications.



### 5 Analysis

CFX Maestro Software provides intuitive setup and data analysis for all your quantitative PCR needs.



\* U.S. patents 6,627,424; 7,541,170; and 7,560,260.

## Ordering Information

Catalog # Description

### Cell Lysis and RNA Isolation

<b>1</b>	7326820	<b>Aurum Total RNA Mini Kit</b> , 50 preps, includes 50 RNA binding columns, 50 capless collection tubes (2.0 ml), 100 capped sample tubes (2.0 ml), 50 capped sample tubes (1.5 ml), 1 vial lyophilized DNase I, RNase-free reagents
	7326830	<b>Aurum Total RNA Fatty and Fibrous Tissue Kit</b> , 50 preps, includes 50 RNA binding columns, 50 capless collection tubes (2.0 ml), 100 capped sample tubes (2.0 ml), 50 capped sample tubes (1.5 ml), 1 vial lyophilized DNase I, RNase-free reagents, plasticware, 50 ml PureZOL RNA Isolation Reagent
	7326800	<b>Aurum Total RNA 96 Kit</b> , 2 x 96-well preps, includes 2 growth membranes, sealing tape, 2 RNA binding plates, 2 collection microplates, 2 vials lyophilized DNase I, RNase-free reagents
	7326880	<b>PureZOL RNA Isolation Reagent</b> , 50 ml, reagent for DNA- and protein-free RNA isolation
	1725085	<b>SingleShot SYBR® Green Kit for Cell Lysis and RT-qPCR</b> , 100 x 50 µl reactions
	1725090	<b>SingleShot Probes Kit for Cell Lysis and RT-qPCR</b> , 100 x 50 µl reactions

More at [bio-rad.com/RNAIsolation](http://bio-rad.com/RNAIsolation)

### Target Selection

#### PrimePCR Primers, Assays, and Arrays

Bio-Rad offers bench-tested primers and probes in predesigned plate, custom plate, and individual assay formats.

More at [bio-rad.com/PrimePCR](http://bio-rad.com/PrimePCR)

### Reaction Setup

<b>3</b>	<b>iScript, SsoAdvanced, Reliance, and iTaq Supermixes</b>	
	1708841	<b>iScript Reverse Transcription Supermix</b> , 400 µl (4 x 100 µl vials), 100 x 20 µl reactions
	1725035	<b>iScript gDNA Clear cDNA Synthesis Kit</b> , 100 x 20 µl reactions
	1725038	<b>iScript Advanced cDNA Synthesis Kit</b> , 100 x 20 µl reactions
	1708891	<b>iScript cDNA Synthesis Kit</b> , 100 x 20 µl reactions
	1725271	<b>SsoAdvanced Universal SYBR® Green Supermix</b> , 5 ml (5 x 1 ml vials), 500 x 20 µl reactions
	1725281	<b>SsoAdvanced Universal Probes Supermix</b> , 5 ml (5 x 1 ml vials), 500 x 20 µl reactions
	1725017	<b>SsoAdvanced Universal Inhibitor-Tolerant SYBR® Green Supermix</b> , 5 ml (1 x 5 ml vial), 500 x 20 µl reactions
	1725160	<b>SsoAdvanced PreAmp Supermix</b> , 1.25 ml (1 x 1.25 ml vial), 50 x 50 µl reactions
	12010220	<b>Reliance One-Step Multiplex RT-qPCR Supermix</b> , 5 ml (5 x 1 ml vials), 1,000 x 20 µl reactions
	12010221	<b>Reliance One-Step Multiplex RT-qPCR Supermix</b> , 10 ml (2 x 5 ml vials), 2,000 x 20 µl reactions
	1725120	<b>iTaq Universal SYBR® Green Supermix</b> , 2 ml (2 x 1 ml vials), 200 x 20 µl reactions
	1725151	<b>iTaq Universal SYBR® Green One-Step Kit</b> , 500 x 20 µl reactions
	1725130	<b>iTaq Universal Probes Supermix</b> , 2 ml (2 x 1 ml vials), 200 x 20 µl reactions
	1725141	<b>iTaq Universal Probes One-Step Kit</b> , 500 x 20 µl reactions

More at [bio-rad.com/PCRReagents](http://bio-rad.com/PCRReagents)

Catalog # Description

### PCR Plastics and Sealer

<b>3</b>	HSP9601	<b>Hard-Shell Low-Profile 96-Well Skirted PCR Plates</b> , pkg of 50, white shell/clear well, rigid 2-component design
	HSP3805	<b>Hard-Shell 384-Well Skirted PCR Plates</b> , pkg of 50, clear shell/white well for higher fluorescent signals, rigid 2-component design
	1814000	<b>PX1 PCR Plate Sealer</b> , includes heat sealing instrument, plate support block that holds 96-well and 384-well plates, sealing frame, power cord
	MSB1001	<b>Microseal 'B' Adhesive Seals</b> , optically clear, 100
	MSF1001	<b>Microseal 'F' PCR Strip Seals</b> , foil, pierceable, 100
	TCS0803	<b>Optical Flat 8-Cap Strips</b> , for 0.2 ml PCR tubes and plates, ultraclear, 120

More at [bio-rad.com/PCRPlastics](http://bio-rad.com/PCRPlastics)

### Detection

<b>4</b>	12011319	<b>CFX Opus 96 Real-Time PCR System</b> , includes integrated 96-well system, Ethernet cable, USB cable, and power cord
	17005940	<b>CFX Opus 96 Real-Time PCR System with Starter Pack</b> , includes integrated 96-well system; Ethernet cable; USB cable; power cord; Hard-Shell 96-Well PCR Plates, low profile, thin wall, skirted, white well/clear shell; SsoAdvanced Universal SYBR® Green Supermix, 2 x 1 ml; Microseal 'B' Adhesive Seals, 100; CFX Maestro Software
	12011452	<b>CFX Opus 384 Real-Time PCR System</b> , includes integrated 384-well system, Ethernet cable, USB cable, and power cord
	17005939	<b>CFX Opus 384 Real-Time PCR System with Starter Pack</b> , includes integrated 384-well system; Ethernet cable; USB cable; power cord; Hard-Shell 384-Well PCR Plates, thin wall, skirted, clear shell/white well, barcoded; SsoAdvanced Universal SYBR® Green Supermix, 2 x 1 ml; Microseal 'B' Adhesive Seals, 100; CFX Maestro Software
	1854095	<b>CFX96 Touch Deep Well Real-Time PCR System</b> , includes C1000 Touch Thermal Cycler Chassis, CFX96 Deep Well Optical Reaction Module, cables
	1855201	<b>CFX Connect Real-Time PCR System</b> , includes CFX Connect Thermal Cycler Chassis, CFX Connect Optical Reaction Module, USB cable

More at [bio-rad.com/qPCR](http://bio-rad.com/qPCR)

### Analysis

	All software for Windows unless otherwise noted.	
	12013758	<b>CFX Maestro Software 2.0</b>
	12004128	<b>CFX Maestro Software for Mac</b>
	12012832	<b>CFX Maestro Software 2.0, Security Edition</b> , 1 license
	12013028	<b>CFX Maestro Software 2.0, Security Edition</b> , 5 licenses
	12012834	<b>CFX Maestro Software 2.0, Chinese Edition</b>
	12012833	<b>CFX Maestro Software 2.0, Russian Edition</b>
	1845025	<b>Precision Melt Analysis Software</b> , 2 user licenses, data analysis software for all CFX Real-Time PCR Systems, 2 HASP HL keys, calibration kit

More at [bio-rad.com/CFXMaestro](http://bio-rad.com/CFXMaestro)

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

All trademarks used herein are the property of their respective owner.

SYBR is a trademark of Thermo Fisher Scientific Inc. Bio-Rad Laboratories, Inc. is licensed by Thermo Fisher Scientific Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Hard-Shell Plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Website [bio-rad.com](http://bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

