

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.

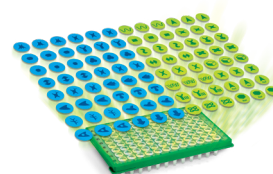
1 Cell Lysis and RNA Isolation

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 — Sso7d fusion polymerase technology provides robust performance.

Plastics — 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.



Ordering Information

Catalog # Description

Cell Lysis and RNA Isolation

7326820	Aurum Total RNA Mini Kit , 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	Aurum Total RNA Fatty and Fibrous Tissue Kit , 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	Aurum Total RNA 96 Kit , 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	PureZOL RNA Isolation Reagent , 50 ml, reagent for DNA- and protein-free RNA isolation
1725085	SingleShot SYBR® Green Kit for Cell Lysis and RT-qPCR , 100 x 50 µl reactions
1725090	SingleShot Probes Kit for Cell Lysis and RT-qPCR , 100 x 50 µl reactions

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

Target Selection

PrimePCR Primers, Assays, and Arrays

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

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

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.
© 2025 Bio-Rad Laboratories, Inc.

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.

Catalog # Description

Reaction Setup

iScript, SsoAdvanced, Reliance, and iTaq Supermixes

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

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

PCR Plastics and Sealer

HSP9601	Hard-Shell Low-Profile 96-Well Skirted PCR Plates , pkg of 50, white shell/clear well, rigid 2-component design
HSP3805	Hard-Shell 384-Well Skirted PCR Plates , pkg of 50, clear shell/white well for higher fluorescent signals, rigid 2-component design
1814000	PX1 PCR Plate Sealer , includes heat sealing instrument, plate support block that holds 96-well and 384-well plates, sealing frame, power cord

Catalog # Description

MSB1001B	Microseal™ 'B' Adhesive Seals , optically clear, 100
MSF1001B	Microseal 'F' PCR Plate Seals , foil, pierceable, 100
TCS0803	Optical Flat 8-Cap Strips , for 0.2 ml PCR tubes and plates, ultraclear, 120

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

Detection

12011319	CFX Opus 96 Real-Time PCR System , includes integrated 96-well system, Ethernet cable, USB cable, and power cord
17005940	CFX Opus 96 Real-Time PCR System with Starter Pack , 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	CFX Opus 384 Real-Time PCR System , includes integrated 384-well system, Ethernet cable, USB cable, and power cord
17005939	CFX Opus 384 Real-Time PCR System with Starter Pack , 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
12016658	CFX Opus Deepwell Real-Time PCR System , includes integrated 96-well sample block for sample reactions of 10–125 µl, ethernet cable, USB cable, and power cord
12016265	CFX Duet Real-Time PCR System , includes fixed 96-well sample block, 2-color multiplex and FRET detection, ethernet cable, USB cable, and power cord

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

Analysis

All software for Windows unless otherwise noted.

12013758	CFX Maestro Software
12012832	CFX Maestro Software, Security Edition , 1 license
12013028	CFX Maestro Software, Security Edition , 5 licenses
12012834	CFX Maestro Software, Chinese Edition
12012833	CFX Maestro Software, Russian Edition

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



Bio-Rad
Laboratories, Inc.

Life Science Group

Website [bio-rad.com](https://www.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 1 800 361 1808 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 82 080 007 7373 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