



## 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

### Isolate RNA

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



2

### Synthesize cDNA

iScript™ Reverse Transcription Kits provide simple, fast, and sensitive first-strand cDNA synthesis.



3

### Set Up the Plate

**PrimePCR™ Assays** — gene-specific primers are bench tested for proven performance.

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

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



4

### Run Real-Time PCR

Bio-Rad's real-time PCR detection systems feature advanced optical technology and precise thermal control for optimal performance.



5

### Analyze Data

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
<b>1</b> 7326820	<b>Isolate RNA</b> <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

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

Catalog #	Description
<b>2</b> 1708840	<b>Synthesize cDNA</b> <b>iScript Reverse Transcription Supermix for RT-qPCR</b> , 25 x 20 µl reactions, 4 x 25 µl 5x supermix contains reverse transcriptase, RNase inhibitor, dNTPs, primers, MgCl <sub>2</sub> , stabilizers; 50 reactions of no-RT control supermix
1725037	<b>iScript Advanced cDNA Synthesis Kit for RT-qPCR</b> , 25 x 20 µl reactions, includes 100 µl 5x iScript Advanced Reaction Mix, 25 µl iScript Advanced Reverse Transcriptase, nuclease-free water
1708890	<b>iScript cDNA Synthesis Kit</b> , 25 x 20 µl reactions, includes 100 µl 5x iScript Reaction Mix, 25 µl iScript Reverse Transcriptase, nuclease-free water

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

Catalog #	Description
<b>3</b>	<b>Set Up the Plate</b> <b>PrimePCR Assays</b> Bio-Rad offers bench-tested primers and probes in predesigned plate, custom plate, and individual assay formats. Go to <a href="http://bio-rad.com/web/PrimePCRAssays">bio-rad.com/web/PrimePCRAssays</a> for more information.
	<b>SsoAdvanced™ and iTaq™ Supermixes</b>
1725270	<b>SsoAdvanced™ Universal SYBR® Green Supermix</b> , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, Sso7d fusion polymerase, MgCl <sub>2</sub> , SYBR® Green I, ROX Normalization Dyes
1725280	<b>SsoAdvanced Universal Probes Supermix</b> , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, Sso7d fusion polymerase, MgCl <sub>2</sub> , stabilizers, ROX Normalization Dyes
1725120	<b>iTaq™ Universal SYBR® Green Supermix</b> , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, iTaq DNA Polymerase, MgCl <sub>2</sub> , SYBR® Green I, enhancers, stabilizers, fluorescein, and ROX Normalization Dyes
1725130	<b>iTaq Universal Probes Supermix</b> , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, iTaq DNA Polymerase, MgCl <sub>2</sub> , enhancers, stabilizers, ROX Normalization Dyes

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

Catalog #	Description
<b>3</b> HSP9601	<b>PCR Plates and Sealer</b> <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

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

Catalog #	Description
<b>4</b> 1855195	<b>Run Real-Time PCR</b> <b>CFX96 Touch™ Real-Time PCR Detection System</b> , includes C1000 Touch™ Thermal Cycler Chassis, CFX96™ Optical Reaction Module, USB cable
1855485	<b>CFX384 Touch™ Real-Time PCR Detection System</b> , includes C1000 Touch Thermal Cycler Chassis, CFX384™ Optical Reaction Module, USB cable
1855201	<b>CFX Connect™ Real-Time PCR Detection System</b> , includes CFX Connect Thermal Cycler Chassis, CFX Connect Optical Reaction Module, USB cable

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

Catalog #	Description
<b>5</b> 12004110	<b>Analyze Data</b> <b>CFX Maestro Software</b> , Windows PC version. Mac, Security Edition (U.S. FDA 21 CFR Part 11 compliant), and Russian and Chinese language versions are also available

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

Mac is a trademark of Apple Inc. ROX and SYBR are trademarks 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. Windows is a trademark of Microsoft Corporation.

Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

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

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

