



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 Manager™ 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

1	Catalog #	Description
1	Isolate RNA 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
More at bio-rad.com/web/RNASolutions		
2	Synthesize cDNA 1708840	iScript Reverse Transcription Supermix for RT-qPCR , 25 x 20 µl reactions, 4 x 25 µl 5x supermix contains reverse transcriptase, RNase inhibitor, dNTPs, primers, MgCl ₂ , stabilizers; 50 reactions of no-RT control supermix
	1725037	iScript Advanced cDNA Synthesis Kit for RT-qPCR , 25 x 20 µl reactions, includes 100 µl 5x iScript Advanced Reaction Mix, 25 µl iScript Advanced Reverse Transcriptase, nuclease-free water
	1708890	iScript cDNA Synthesis Kit , 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		
3	Set Up the Plate PrimePCR Assays	Bio-Rad offers bench-tested primers and probes in predesigned plate, custom plate, and individual assay formats. Go to bio-rad.com/web/PrimePCRAssays for more information.
	SsoAdvanced™ and iTaq™ Supermixes	
	1725270	SsoAdvanced™ Universal SYBR® Green Supermix , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , SYBR® Green I, ROX Normalization Dyes
	1725280	SsoAdvanced Universal Probes Supermix , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , stabilizers, ROX Normalization Dyes
	1725120	iTaq™ Universal SYBR® Green Supermix , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, iTaq DNA Polymerase, MgCl ₂ , SYBR® Green I, enhancers, stabilizers, fluorescein, and ROX Normalization Dyes
1725130	iTaq Universal Probes Supermix , 200 x 20 µl reactions, 2 ml (2 x 1 ml vials), 2x qPCR mix, contains dNTPs, iTaq DNA Polymerase, MgCl ₂ , enhancers, stabilizers, ROX Normalization Dyes	
More at bio-rad.com/web/PCRSupermixes		

3	Catalog #	Description
3	PCR Plates 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
More at bio-rad.com/web/PlasticSolutions		
4	Run Real-Time PCR 1855195	CFX96 Touch™ Real-Time PCR Detection System , includes C1000 Touch™ Thermal Cycler Chassis, CFX96™ Optical Reaction Module, CFX Manager Software, license for qbase+ Software, communication cable
	1855485	CFX384 Touch™ Real-Time PCR Detection System , includes C1000 Touch Thermal Cycler Chassis, CFX384™ Optical Reaction Module, CFX Manager Software, license for qbase+ Software, communication cable
	1855201	CFX Connect™ Real-Time PCR Detection System , includes CFX Connect Thermal Cycler Chassis, CFX Connect Optical Reaction Module, CFX Manager Software, license for qbase+ Software, communication cable
More at bio-rad.com/web/RealTimePCR		
5	Analyze Data	CFX Manager Software can be obtained with the purchase of any Bio-Rad real-time PCR detection system.
	More at bio-rad.com/web/CFXManager	

ROX and SYBR are trademarks of Life Technologies Corporation. Bio-Rad Laboratories, Inc. is licensed by Life Technologies Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

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.

The use of iTaq and SsoAdvanced Supermixes is covered by one or more of the following U.S. patents and corresponding patent claims outside the U.S.: 5,804,375; 5,994,056; and 6,171,785. The purchase of these products includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. These products are for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

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 www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000