

Gene Expression



Description

Purification of RNA from tissues, cell cultures, or bacterial cells

Analysis of RNA for quality and quantity

Applications

Nucleic acid profiles

Technologies

Nucleic acid sample preparation

Automated nucleic acid electrophoresis
Nucleic acid electrophoresis
Image analysis
Spectrophotometry and fluorometry

Related Products



Aurum™ total RNA kits
PureZOL™ RNA isolation reagent



Experion™ automated electrophoresis system
Experion RNA analysis kits
Wide mini ReadySub-Cell™ GT cell
ReadyAgarose™ gels
Molecular Imager® Gel Doc™ XR system
Quantity One® 1-D analysis software
SmartSpec™ Plus spectrophotometer

Amplification

Reverse transcription and amplification of transcripts

Reverse transcription PCR (RT-PCR)



iQ™5 multicolor real-time PCR detection system

iQ™ supermix

iCycler® thermal cycler

MyCycler™ thermal cycler

iScript™ cDNA synthesis kits

Profiling and Quantitation

Profiling and quantitation of transcripts

Quantitative RT-PCR (RT-qPCR)

Microarray analysis

Nucleic acid blotting

Nucleic acid sequencing

Multiplex suspension array technology — nucleic acid

Image analysis

Data analysis and bioinformatics



MiniOpticon™ two-color real-time PCR detection system

MyiQ™ single-color real-time PCR detection system

BioOdyssey™ Calligrapher™ miniarrayer

Bio-Plex® 200 system with HTF

iTaq™ SYBR® Green supermix with ROX

iQ™ SYBR® Green supermix

iTaq™ DNA polymerase

Molecular Imager Gel Doc XR system

Quantity One 1-D analysis software


Gene Expression Profile

From RNA purification to profiling and quantitation, Bio-Rad provides life science researchers with superior tools, service, and support that produce consistently reliable and sensitive results for characterizing gene expression.

Explore the transcriptome with Bio-Rad's full line of genomic reagents and instruments. Whether you are analyzing thousands of genes or just one, Bio-Rad Laboratories has the equipment to meet your needs. Learn how our reagents and instrumentation can transform your laboratory and facilitate your workflow. Visit us on the Web at www.bio-rad.com/genomics/

For a complete list of products and related product information, refer to the 2006/2007 catalog, visit us on the Web at www.bio-rad.com, or contact your local Bio-Rad representative.

Ordering Information

Catalog #	Description	Catalog #	Description
Sample Preparation			
732-6800	Aurum Total RNA 96 Kit, 2 x 96-well preps	169-2000	BioOdyssey Calligrapher MiniArrayer, 115 V
732-6890	PureZOL RNA Isolation Reagent, 100 ml	169-2100	BioOdyssey Calligrapher MiniArrayer, 230 V
		169-2200	BioOdyssey Calligrapher MiniArrayer With Cooling Module, 115 V
Sample Assessment			
700-7001	Experion System, for RNA analysis	169-2300	BioOdyssey Calligrapher MiniArrayer With Cooling Module, 230 V
700-7103	Experion RNA StdSens Analysis Kit	165-5003	Model 185 Vacuum Blotter Basic Unit
700-7105	Experion RNA HighSens Analysis Kit	Varies	Sequi-Gen® GT Systems
170-4489	Wide Mini ReadySub-Cell GT Cell	170-9300	FPQuest™ Software
161-3049	Wide Mini ReadyAgarose Gel, TAE, 1.0%, 32-well	171-000201	Bio-Plex 200 System
161-3053	Wide Mini ReadyAgarose Gel, TAE, 1.0%, 2 x 32-well	171-000205	Bio-Plex 200 System With HTF, includes high-throughput fluidics (HTF)
161-3050	Wide Mini ReadyAgarose Gel, TAE, 1.0% plus ethidium bromide, 32-well	Varies	Bio-Plex Assays
161-3054	Wide Mini ReadyAgarose Gel, TAE, 1.0% plus ethidium bromide, 2 x 32-well	170-8850	iTaq SYBR Green Supremix With ROX, 200 x 50 µl reactions, 2x mix contains 0.4 mM each of dATP, dCTP, and dGTP, 0.8 mM dUTP, 50 U/ml iTaq DNA polymerase, 6 mM Mg ²⁺ , SYBR Green I, ROX reference dye, stabilizers
170-8170	Molecular Imager Gel Doc XR System, PC, includes Quantity One software	170-8880	iQ SYBR Green Supremix, 100 x 50 µl reactions
170-8171	Molecular Imager Gel Doc XR System, Mac, includes Quantity One software	170-8870	iTaq DNA Polymerase, 5 U/µl, includes 250 U polymerase, 1.25 ml of 10x PCR buffer, 1.25 ml of 50 mM MgCl ₂ solution
170-9600	Quantity One 1-D Analysis Software		
170-2525	SmartSpec Plus Spectrophotometer		
170-2402	VersaFluor Fluorometer		
Amplification			
170-9780	iQ5 Multicolor Real-Time PCR Detection System, includes iCycler base, optics module, software CD-ROM, 5 installed filter sets, 96-well reaction module, calibration solutions, optical-quality 96-well PCR plates, Microseal 'B' seals, communication cables, power cord, quick reference cards, instructions		LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only.
170-8860	iQ Supremix, 100 x 50 µl reactions, 2x mix		Certain arrays and/or the method of preparation, analysis, or use may be covered by patents or other intellectual property rights owned by third parties.
170-8720	iCycler Thermal Cycler With 96 x 0.2 ml Reaction Module, includes iCycler base with 96 x 0.2 ml iCycler reaction module, in-sample temperature probe (0.2 ml tube size), PCR tubes, power cord, instructions, quick-reference card		The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.
170-9703	MyCycler Thermal Cycler System, includes MyCycler thermal cycler, PCR tubes, iTaq DNA polymerase, power cord, quick reference guide, instructions		SYBR is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.
170-8894	iScript One-Step RT-PCR Kit for Probes, 50 x 50 µl reactions		Notice regarding Bio-Rad thermal cyclers and real-time systems: Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents Nos. 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160-163 and 167 only), and 6,703,236 (claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applera Corporation. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.
170-8892	iScript™ One-Step RT-PCR Kit With SYBR® Green, 50 x 50 µl reactions		Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics.
Profiling and Quantitation			
CFB-3120	MiniOpticon Real-Time PCR System, includes optical housing, MJ Mini™ thermal cycler, analysis software		
CFD-3121	MiniOpticon Real-Time PCR Detector, for MJ Mini cycler upgrade		
170-9770	MiQ Single-Color Real-Time PCR Detection System, includes iCycler base, optics module, software CD-ROM, 96-well optical reaction module, optical-quality 96-well PCR plates, Microseal 'B' seals, communication cables, power cords, instructions		



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500 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 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000