



Life Science Research  
Biomarker Discovery



# Biomarker Discovery





**Bioinformatics**

**Assay Development**

**Validation**

**Assay for Disease State**

Analysis of profiling and expression data

Development of assays for biomarker(s) for disease states

Validation of specific biomarker assays

From sample preparation to validation, Bio-Rad's uniquely comprehensive suite of tools, service, and support deliver consistent, reliable results to drive the biomarker discovery efforts of life science researchers.

Bio-Rad's genomics and proteomics tools offer complete application solutions, including innovative products that facilitate the biomarker discovery workflow. For more information on how Bio-Rad can help you achieve your research goals, visit us on the Web at [www.bio-rad.com](http://www.bio-rad.com)

Data analysis and bioinformatics

**Protein**

**Nucleic Acid**

**Protein**

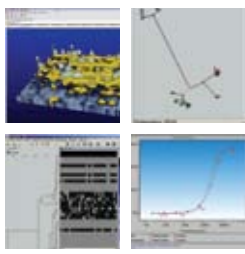
**Nucleic Acid**

- Protein blotting
- Multiplex suspension array technology — protein
- Surface-enhanced laser desorption/ionization (SELDI)
- ELISA
- Spectrometry and fluorometry

- Nucleic acid blotting
- Quantitative PCR (qPCR)

- Protein interaction analysis
- Multiplex suspension array technology — protein
- Microarray analysis
- Protein blotting

- Microarray analysis
- Nucleic acid blotting
- Quantitative PCR (qPCR)



- Proteomweaver™ 2-D analysis software
- FPQuest™ software
- InfoQuest™ FP software
- Bio-Plex Manager™ 4.1 Standard Edition software




- Trans-Blot® Plus cell
- Bio-Plex 200 system
- Ultramark™ microplate system
- ProteinChip SELDI system
- iQ™5 multicolor real-time PCR detection system



- ProteOn™ XPR36 protein interaction array system
- BioOdyssey Calligrapher miniarrayer
- MiniOpticon two-color real-time PCR detection system
- Criterion blotter
- PowerPac HC power supply

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](http://www.bio-rad.com), or contact your local Bio-Rad representative.

**Ordering Information**

Catalog #	Description	Catalog #	Description
<b>Sample Preparation</b>			
732-6800	<b>Aurum Total RNA 96 Kit</b> , 2 x 96-well preps	169-2000	<b>BioOdyssey Calligrapher MiniArrayer</b> , 115 V
163-2086	<b>ReadyPrep Protein Extraction Kit (Total Protein)</b> , 20 preps	169-2100	<b>BioOdyssey Calligrapher MiniArrayer</b> , 230 V
163-2100	<b>ReadyPrep Sequential Extraction Kit</b> , 5–15 preps	CFB-3120	<b>MiniOpticon Real-Time PCR System</b> , includes optical housing, MJ Mini thermal cycler, analysis software
163-2130	<b>ReadyPrep 2-D Cleanup Kit</b> , 50 preps	CFD-3121	<b>MiniOpticon Real-Time PCR Detector</b> , for MJ Mini cycler upgrade
732-6340	<b>AquaPure™ Genomic DNA Isolation Kit</b> , for cultured cells and gram-negative bacteria	<b>Bioinformatics</b>	
142-2825	<b>Chelex® 100 Resin</b> , 100 g, iron, 100–200 dry mesh, 150–300 µm wet bead	165-9770	<b>Proteomweaver 4.0 Professional 2-D Analysis Software</b>
<b>Sample Assessment</b>			
165-6001	<b>Criterion Cell</b>	170-9300	<b>FPQuest Software</b>
Varies	<b>Criterion Precast Gels</b>	170-9310	<b>InfoQuestFP Basic Fingerprint Types Module</b>
700-7000	<b>Experion System</b> , for protein analysis (analysis kits sold separately)	170-9311	<b>InfoQuestFP Character Types Module</b>
700-7001	<b>Experion System</b> , 100–120 V, for RNA analysis (analysis kits sold separately)	170-9312	<b>InfoQuestFP Sequence Types Module</b>
700-7002	<b>Experion System</b> , 220–240 V, for RNA analysis (analysis kits sold separately)	170-9313	<b>InfoQuestFP Matrix Types Module</b>
700-7102	<b>Experion Pro260 Analysis Kit for 25 Chips</b>	171-STND01	<b>Bio-Plex Manager 4.1 Standard Software</b> , 1-user desktop license
700-7104	<b>Experion RNA StdSens Analysis Kit for 25 Chips</b> , includes 25 RNA StdSens chips, Experion RNA StdSens reagents and supplies for 25 chips	171-SCRT00	<b>Bio-Plex Manager 4.1 Security Workstation License</b>
700-7106	<b>Experion RNA HighSens Analysis Kit for 25 Chips</b> , includes 25 RNA HighSens chips, Experion RNA HighSens reagents and supplies for 25 chips	<b>Assay Development</b>	
165-4150	<b>PROTEAN Plus Dodeca Cell</b> , 100/120 V	170-3990	<b>Trans-Blot Plus Cell With Plate Electrodes and Super Cooling Coil</b>
165-4151	<b>PROTEAN Plus Dodeca Cell</b> , 220/240 V	170-9500	<b>Ultramark Microplate Imaging System</b>
164-5050	<b>PowerPac Basic Power Supply</b>	170-9549	<b>Ultramark™ EX Microplate Imaging System</b>
161-0491	<b>Flamingo Fluorescent Gel Stain</b> , 10x solution, 100 ml	170-9780	<b>iQ5 Multicolor Real-Time PCR Detection System</b> , 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
170-2525	<b>SpartSpec Plus Spectrophotometer</b>	<b>Validation</b>	
170-2402	<b>VersaFluor Fluorometer</b>	176-0100	<b>ProteOn XPR36 Protein Interaction Array System</b>
500-0205	<b>Quick Start™ Bradford 1x Dye Reagent</b> , 1 L	170-4070	<b>Criterion Blotter With Plate Electrodes</b>
<b>Profiling and Quantitation</b>			
171-000201	<b>Bio-Plex 200 System</b>	164-5052	<b>PowerPac HC Power Supply</b>
171-000205	<b>Bio-Plex 200 System With HTF</b> , includes high-throughput fluidics (HTF)	 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.	
Z50-00011	<b>ProteinChip SELDI System</b> , Enterprise Edition, 115 V	The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.	
Z50-00013	<b>ProteinChip SELDI System</b> , Personal Edition, 115 V	The SELDI process is covered by U.S. Patents 5,719,060, 6,225,047, 6,579,719, 6,818,411 and other issued patents and pending applications in the U.S. and other jurisdictions.	
Z50-00021	<b>ProteinChip SELDI System</b> , Enterprise Edition, 230 V	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.	
Z50-00023	<b>ProteinChip SELDI System</b> , Personal Edition, 230 V	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.	
Varies	<b>ProteinChip Arrays</b>		
Varies	<b>ProteinChip Energy Absorbing Molecules (EAMs)</b>		
C10-00001	<b>ProteinChip Protein Calibrant Kit</b>		
C10-00002	<b>ProteinChip Peptide Calibrant Kit</b>		
170-8640	<b>Molecular Imager VersaDoc MP 4000 System</b> , includes Quantity One software		
170-9460	<b>Molecular Imager PharosFX Plus System</b> , includes Quantity One software		
165-7200	<b>EXQuest Spot Cutter</b>		



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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