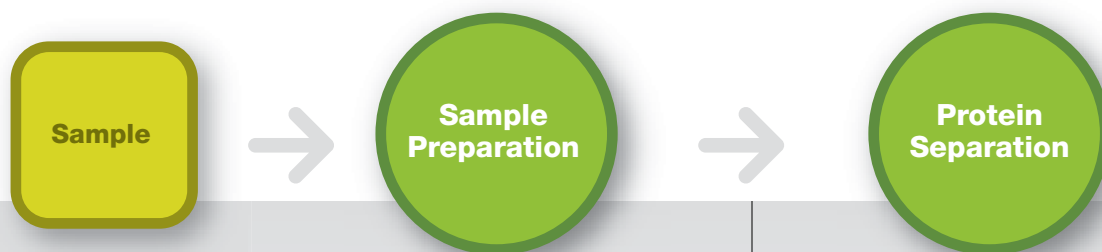




Life Science Research
Protein Separation and Analysis



Protein Separation and Analysis



Description

Protein purification and fractionation from tissues, cell cultures, or bacterial cells

Separation of proteins by electrophoresis, chromatography, or a combination of methods

Applications

- Ion exchange for protein identification or profiling
- Affinity and 1-D electrophoresis for protein identification or profiling
- Chromatography and electrophoresis for protein identification or profiling
- 1-D electrophoresis for detection or quantitation
- 2-D electrophoresis for detection or quantitation

-
-
-
-
-

Technologies

Protein sample preparation

Chromatography

- Ion exchange chromatography
- Hydroxyapatite and affinity chromatography
- Size exclusion chromatography
- Automated affinity chromatography

Electrophoresis

- Protein electrophoresis
- 2-D electrophoresis
- Isoelectric focusing
- Preparative electrophoresis

Related Products



- ReadyPrep™ protein extraction kits
- ReadyPrep sequential extraction kit
- Model 491 prep cell
- MicroRotor™ cell



- BioLogic DuoFlow™ F10 workstation
- Profinia™ purification system
- UNOsphere™ Q support
- CHT™ ceramic hydroxyapatite
- Bio-Gel® P polyacrylamide gels
- PROTEAN® IEF system
- Mini-PROTEAN® Tetra cell
- Criterion™ cell and precast gels
- PROTEAN® II XL cell
- PowerPac™ Basic power supply



Protein profiling, characterization, quantitation, and/or identification

Analyzed Proteins

For researchers who are discovering and characterizing proteins, Bio-Rad's protein analysis tools offer a complete system approach, a unique position among technology suppliers, that delivers high-value, accessible technologies.

From sample preparation to protein analysis and profiling, Bio-Rad's protein separation and analysis workflow tools provide choices in methodology, protocols, and products. The protein separation and analysis workflow focuses on five technologies: sample preparation, chromatography, electrophoresis, imaging and analysis, and surface-enhanced laser desorption/ionization (SELDI). Each technology is well suited to the separation, analysis, and profiling of protein samples.

Identification

- Gel excision
- Matrix-assisted laser desorption/ionization (MALDI)
- Electrospray ionization (ESI)

Profiling

- Surface-enhanced laser desorption/ionization (SELDI)
- Multiplex suspension array technology — protein
- Microarray analysis

Characterization and Quantitation

- Image analysis
- Protein blotting
- Microarray analysis
- Data analysis and bioinformatics
- Multiplex suspension array technology — protein
- Protein assays



- EXQuest™ spot cutter
- ProteinChip® SELDI system
- Bio-Plex® 200 system with HTF
- BioOdyssey™ Calligrapher™ miniarrayer
- Trans-Blot® Plus cell
- Molecular Imager® VersaDoc™ MP 4000 system
- Molecular Imager® PharosFX™ system

Protein Separation and Analysis

Sample > Sample Preparation > Protein Separation > Profiling and Quantitation > Analyzed Proteins

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			
163-2086	ReadyPrep Protein Extraction Kit (Total Protein), 20 preps	Varies	Ready Gel Precast Gels
163-2085	ReadyPrep Protein Extraction Kit (Soluble/ Insoluble), 20 preps	165-6001	Criterion Cell
163-2089	ReadyPrep Protein Extraction Kit (Cytoplasmic/ Nuclear), 50 preps	Varies	Criterion Precast Gels
163-2088	ReadyPrep Protein Extraction Kit (Membrane I), 50 preps	165-3188	PROTEAN II XL Cell, wide-format
163-2130	ReadyPrep 2-D Cleanup Kit, 50 preps	161-0375	Precision Plus Protein Kaleidoscope Standards, 500 µl, 50 applications
163-2100	ReadyPrep Sequential Extraction Kit, 5–15 preps	164-5050	PowerPac Basic Power Supply
170-2926	Model 491 Prep Cell, 100/120 V	Profiling and Quantitation	
170-2927	Model 491 Prep Cell, 220/240 V	165-7200	EXQuest Spot Cutter
170-2800	MicroRotor Cell Kit, 100/120 V	165-7201	EXQuest Spot Cutter With PC
170-2801	MicroRotor Cell Kit, 220/240 V	Z50-00011	ProteinChip SELDI System, Enterprise Edition, 115 V
170-2802	MicroRotor System, 100/120 V, includes kit with PowerPac HV power supply	Z50-00013	ProteinChip SELDI System, Personal Edition, 115 V
170-2803	MicroRotor System, 220/240 V	Z50-00021	ProteinChip SELDI System, Enterprise Edition, 230 V
Protein Separation			
760-0037	BioLogic DuoFlow 10 System, 100/120 V, includes Dell controller and monitor, USB Bitbus communicator, F10 workstation, MX-1 mixer, 3-tray rack, AVR7-3 sample inject valve, fittings kit, UV detector with 5 mm flow cell and 254/280 nm filters, conductivity monitor, starter kit, UNO Q1 column, instructions	Z50-00023	ProteinChip SELDI System, Personal Edition, 230 V
760-0038	BioLogic DuoFlow 10 System, 220/240 V	Varies	ProteinChip Arrays
620-1004	Profinia Instrument With Accessory Kit, includes cleaning tray, inline filter pack, 2 x 50 ml sample lids, 2 x 15 ml sample lids, bottle starter pack, waste/diluent bottle set	Varies	ProteinChip Energy Absorbing Molecules (EAMs)
620-1005	Profinia Instrument With Accessory Kit and Native IMAC Starter Kit, includes same as 620-1004 with Profinia native IMAC buffer kit, 1 x 1 ml IMAC and 1 x 10 ml desalting cartridge, <i>E. coli</i> lysate	C10-00001	ProteinChip Protein Calibrant Kit
620-1006	Profinia Instrument With Accessory Kit and GST Starter Kit, includes same as 620-1004 with Profinia GST buffer kit, 1 x 1 ml GST and 1 x 10 ml desalting cartridge, <i>E. coli</i> lysate, glutathione reagent	C10-00002	ProteinChip Peptide Calibrant Kit
620-1009	Profinia Protein Purification System, includes same as 620-1004 with Profinia software	171-000201	Bio-Plex 200 System
620-1010	Profinia Protein Purification System With Native IMAC Starter Kit, includes same as 620-1005 with Profinia software	171-000205	Bio-Plex 200 System With HTF, includes high-throughput fluidics (HTF)
620-1011	Profinia Protein Purification System With GST Starter Kit, includes same as 620-1006 with Profinia software	Varies	Bio-Plex Assays
156-0105	UNOsphere Q Support, 500 ml	169-2000	BioOdyssey Calligrapher MiniArrayer, 115 V
158-2000	CHT Ceramic Hydroxyapatite, Type I, 20 µm, 10 g	170-3990	Trans-Blot Plus Cell With Plate Electrodes and Super Cooling Coil
Varies	Bio-Gel P Polyacrylamide Gels	170-8640	Molecular Imager VersaDoc MP 4000 System, PC or Mac, includes Quantity One software
165-4000	PROTEAN IEF System, 90–240 VAC	170-9460	Molecular Imager PharosFX Plus System, PC or Mac, includes Quantity One software
165-4001	PROTEAN IEF Cell, 90–240 VAC, basic unit, includes cell, instructions	170-9450	Molecular Imager PharosFX System, PC or Mac, includes Quantity One software
165-8000	Mini-PROTEAN Tetra Cell, 10-well, 0.75 mm thickness; 4-gel system	165-9770	Proteomweaver 4.0 Professional 2-D Analysis Software
		165-9775	Proteomweaver 4.0 Professional 2-D Analysis Software, with multifluorescence analysis (MFA)
		165-9780	Proteomweaver 4.0 Enterprise 2-D Analysis Software
		165-9790	Proteomweaver 4.0 Enterprise 2-D Analysis Software, with multifluorescence analysis (MFA)
		170-9620	PDQuest Basic 2-D Analysis Software

Mac is a trademark of Apple Corporation.

Purification of fusion proteins may require a license from third parties.

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

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.

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.

BIO-RAD

**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