



Life Science Research
Protein Purification



Protein Purification



Description

Analysis of crude protein extract

Lysis, solubilization, clarification, and fractionation of sample

Analysis of refined protein sample

Applications

- Purification from native secreted protein
- Purification from native intracellular protein
- Purification from recombinant secreted protein
- Purification from recombinant intracellular protein



Technologies

- Automated protein electrophoresis
- Protein electrophoresis
- Protein blotting
- Image analysis
- Spectrophotometry and fluorometry
- Protein assays

Protein sample preparation

- Automated protein electrophoresis
- Protein electrophoresis
- Protein blotting
- Image analysis
- Spectrophotometry and fluorometry
- Protein assays

Related Products



- Experion™ automated electrophoresis system
- Mini-PROTEAN® Tetra cell
- Criterion™ cell and precast gels
- Flamingo™ fluorescent gel stain
- Molecular Imager® Gel Doc™ XR system
- Molecular Imager® ChemiDoc™ XRS system
- SmartSpec™ Plus spectrophotometer
- RC DC™ protein assay



- ReadyPrep™ protein extraction kits
- ReadyPrep sequential extraction kit
- MicroRotofor™ cell



- Flamingo fluorescent gel stain
- SmartSpec Plus spectrophotometer
- RC DC protein assay
- Molecular Imager Gel Doc XR system



Purification of a specific protein or protein family of interest

Assessment of purified protein(s) for quality and quantity

For over half a century, Bio-Rad has provided life science researchers performing protein purification and separation a complete set of easy-to-use, flexible, modular tools that deliver efficient, convenient, and reproducible results.

Bio-Rad offers a complete line of laboratory-scale chromatography instruments, columns, and media that are flexible, upgradable, and easy to use. From automated purification to reliable low-pressure systems, Bio-Rad provides a solution for each phase of purification — capture, intermediate, and polishing — and manufactures media for any scale, from nanograms to kilograms. Visit us on the Web at www.bio-rad.com/chromsystems/

Recombinant

Automated affinity chromatography

Tagged protein purification

Native

Preparative electrophoresis

Ion exchange chromatography

Hydroxyapatite and affinity chromatography

Size exclusion chromatography

Automated protein electrophoresis
 Protein electrophoresis
 Protein blotting
 Image analysis
 Spectrophotometry and fluorometry
 Protein assays



- Profinia™ protein purification system
- Profinity™ IMAC and GST resins
- UNOsphere™ Q support
- Chelex® 100 resin
- Bio-Gel® HT and HTP hydroxyapatite
- CHT™ ceramic hydroxyapatite
- Econo-Pac® protein A columns
- Econo-Column® columns
- BioLogic™ LP system
- Model 491 prep cell

- Experion automated electrophoresis system
- Experion Pro260 analysis kit
- Mini-PROTEAN Tetra cell and Ready Gel® precast gels
- Criterion blotter
- PowerPac™ Universal power supply
- SmartSpec Plus spectrophotometer

Protein Purification

Sample > Sample Assessment > Sample Preparation > Sample Assessment > Purification > Sample Assessment > Purified Protein

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 Assessment		Purification	
700-7060	Experion System With Dell Computer and Monitor , for protein analysis (analysis kits sold separately)	620-1009	Profinia Protein Purification System , 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, with Profinia software
700-7102	Experion Pro260 Analysis Kit for 25 Chips , includes 25 Pro260 chips, Experion Pro260 reagents and supplies for 25 chips	156-0121	Profinity IMAC Uncharged Resin , 10 ml
165-8000	Mini-PROTEANTetra Cell , 10-well, 0.75 mm thickness; 4-gel system	156-0131	Profinity IMAC Ni-Charged Resin , 10 ml
Varies	Ready Gel Precast Gels	732-4612	Bio-Scale Mini Profinity IMAC Cartridge , 1 x 5 ml
165-6001	CriterionCell	732-4622	Bio-Scale Mini Profinity GST Cartridge , 1 x 5 ml
Varies	CriterionTris-HCl Gels	156-0105	UNOsphere Q Support , 500 ml
Varies	CriterionTris-Tricine Gels	142-2825	Chelex 100 Resin , 100 g
165-6024	Criterion Cell/Plate Blotter System , includes Criterion cell and Criterion blotter with plate electrodes	130-0150	Bio-Gel HT Hydroxyapatite , hydrated, 250 ml
164-5070	PowerPac Universal Power Supply	130-0420	Bio-Gel HTP Hydroxyapatite , powder, 100 g
161-0491	Flamingo Fluorescent Gel Stain , 10x solution, 100 ml	130-0520	Bio-Gel HTP Hydroxyapatite , DNA grade, 100 g
161-0787	Bio-Safe™ Coomassie Stain , 5 L	158-2000	CHT Ceramic Hydroxyapatite , 20 µm, Type I, 10 g
170-8170	MolecularImager Gel Doc XR System , PC, includes Quantity One software	158-2200	CHT Ceramic Hydroxyapatite , 20 µm, Type II, 10 g
170-8171	MolecularImager Gel Doc XR System , Mac, includes Quantity One software	Varies	Bio-Gel P Polyacrylamide Gels
170-8650	Molecular Imager® VersaDoc™ MP 5000 System , PC or Mac, includes Quantity One software	732-2022	Econo-PacProtein A Columns , prefilled with Affi-Gel protein A gel, 5
170-8070	Molecular Imager ChemiDoc XRS System , PC, includes Quantity One software	732-1010	Econo-PacChromatography Columns , empty, 50
170-8071	Molecular Imager ChemiDoc XRS System , Mac, includes Quantity One software	732-1011	Econo-PacChromatography Columns , empty, 500
170-2525	SmartSpec Plus Spectrophotometer	Varies	Econo-Column Chromatography Columns
170-2402	VersaFluor Fluorometer	760-0038	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
500-0121	RC DC Protein Assay Kit I , includes RC reagents package, DC protein assay reagents package, bovine γ-globulin standard; 450 standard assays	731-8300	BioLogic DuoFlow 10 System , 220/240 V
Sample Preparation			StandardBioLogic LP System , 100/120 V, includes BioLogic LP controller, BioLogic rack, accessory kit with inject valve (MV-6), proportioning valve/mixer, UV optics, conductivity cell, tubing and fittings kit, column and conductivity cell holder, starter kit, instructions
163-2086	ReadyPrep Protein Extraction Kit (Total Protein) , 20 preps	731-8301	Standard BioLogic LP System , 220/240 V
163-2089	ReadyPrep Protein Extraction Kit (Cytoplasmic/Nuclear) , 50 preps	170-2926	Model491 Prep Cell , 100/120 V
163-2100	ReadyPrepSequential Extraction Kit , 5–15 preps	170-2927	Model491 Prep Cell , 220/240 V
163-2141	MicroRotorforCell Lysis Kit (Mammal) , 15 preps	170-2908	Mini Prep Cell Without Reagent Starter Kit
163-2142	MicroRotorforCell Lysis Kit (Plant) , 10 preps	170-2915	Mini Prep Cell With Reagent Starter Kit
163-2143	MicroRotorforCell Lysis Kit (Yeast) , 15 preps		
163-2144	MicroRotorforCell Lysis Kit (Bacteria) , 15 preps		
170-2800	MicroRotorforCell Kit , 100/120 V		
170-2801	MicroRotorfor Cell Kit , 220/240 V		
170-2802	MicroRotorforSystem , 100/120 V, includes kit with PowerPac HV power supply		
170-2803	MicroRotorforSystem , 220/240 V, includes kit with PowerPac HV power supply		

Coomassie is a trademark of BASF Aktiengesellschaft.



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.

Purification of fusion proteins may require a license from 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