

PROTEIN PURIFICATION Bio-Scale™ Mini Cartridges

- Convenient, prepacked, and disposable design
- Easy connection to liquid chromatography systems using luer lock fittings
- Innovative cartridge design for run-to-run reproducibility
- Wide range of available chromatography media



Convenient, Reliable, and Ready to Use

Chromatography System Compatibility

Bio-Scale Mini cartridges conveniently run a wide variety of chromatography media in the Profinia™ protein purification system, the BioLogic™ LP and BioLogic DuoFlow™ systems, a peristaltic pump, and other chromatography systems. For simple step elution, the cartridges can be used with a luer lock syringe. Fittings are available for connection to HPLC- and FPLC-type chromatography systems.

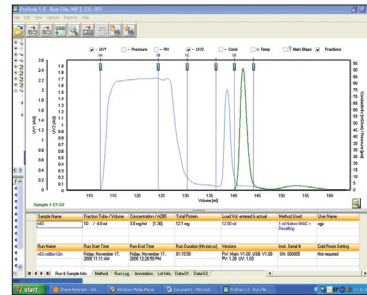
Innovative and Robust Design

Bio-Scale Mini cartridges have a double-wall cartridge design that provides extra durability and allows easy, reliable runs at pressures up to 45 psi and with aqueous buffers most commonly used for protein separation. These new cartridges are robust and constructed of chemically compatible polypropylene. The cartridge design and packing process are a unique patented* Bio-Rad design.

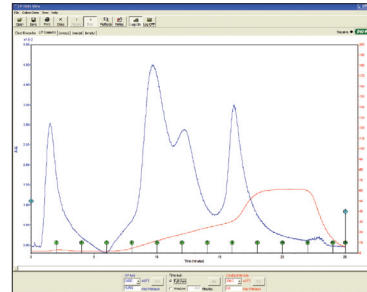
Media Variety and Cartridge Size

The cartridges are prepacked with Bio-Rad's chromatography media for ion exchange, affinity, and size exclusion chromatography. The cartridges come in 1 and 5 ml formats that can be used for method scouting or simple purifications. Larger 10 and 50 ml desalting cartridges are also available. Bio-Scale Mini cartridges are convenient, disposable, and ready to use.

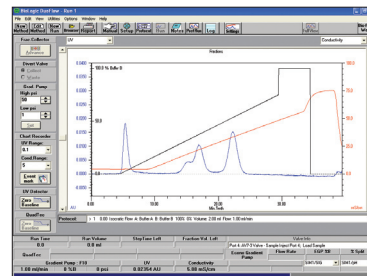
* U.S. patent 7,208,087 B2.



Chromatogram from the Profinia protein purification system using a Bio-Scale™ Mini Profinity™ IMAC cartridge.



Chromatogram from the BioLogic LP system using a Bio-Scale™ Mini Macro-Prep® High Q cartridge.



Chromatogram from the BioLogic DuoFlow system using a Bio-Scale Mini Macro-Prep High S cartridge.



Prepacked Cartridge Selection

Cartridge by Application	Media	Mode of Separation	Chemical Form	Protein Capacity/ml*	Catalog Numbers by Cartridge Size				
					5 x 1 ml	1 x 5 ml	5 x 5 ml		
Protein and Plasmid Purification									
	Macro-Prep High Q	Strong anion exchange	-N ⁺ (CH ₃) ₃	≥37 mg BSA	732-4120	732-4122	732-4124		
	UNOsphere™ Q	Strong anion exchange	-N ⁺ (CH ₃) ₃	180 mg BSA	732-4100	732-4102	732-4104		
	Macro-Prep DEAE	Weak anion exchange	-HN ⁺ (C ₂ H ₅) ₂	>35 mg BSA	732-4140	732-4142	732-4144		
Protein Purification									
	Macro-Prep High S	Strong cation exchange	-SO ₃ ⁻	≥49 mg human IgG	732-4130	732-4132	732-4134		
	UNOsphere Rapid S	Strong cation exchange	-SO ₃ ⁻	60 mg/ml human IgG	732-4400	732-4401	732-4402		
	UNOsphere S	Strong cation exchange	-SO ₃ ⁻	60 mg/ml human IgG	732-4110	732-4112	732-4114		
	Nuvia™ S	Strong cation exchange	-SO ₃ ⁻	110 mg/ml IgG	732-4421	732-4422	732-4423		
Antibody Purification									
	Affi-Gel® Blue DEAE	Affinity	DEAE-Cibacron Blue	0.2–1.0 ml serum F3GA and DEAE	—	732-4632	732-4634		
	UNOsphere SUPra™	Affinity	rProtein A	25–30 mg/ml IgG	732-4201	732-4202	—		
	Affi-Prep® protein A	Affinity	Protein A	8–10 mg mouse monoclonal IgG, 16–23 mg human IgG	732-4600	732-4602	—		
	CHT™ ceramic hydroxyapatite Type I, 40 μm	Mixed mode hydroxyapatite	(Ca ₅ (PO ₄) ₃ OH) ₂	25–60 mg human IgG	—	732-4322	732-4324		
	CHT ceramic hydroxyapatite Type II, 40 μm	Mixed mode hydroxyapatite	(Ca ₅ (PO ₄) ₃ OH) ₂	15–25 mg human IgG	—	732-4332	732-4334		
	CFT™ Ceramic fluoroapatite Type II	Mixed mode fluorapatite	Ca ₁₀ (PO ₄) ₆ F ₂	Lysozyme: 14–21.5 mg/g	—	732-4405	732-4406		
Affinity-Tagged Protein Purification									
	Profinity IMAC	Affinity	Iminodiacetic acid	≥15 mg recombinant histidine (His)-tagged protein	732-4610	732-4612	732-4614		
	Profinity GST	Affinity	Glutathione	≥7 mg/ml recombinant GST-tagged protein	732-4620	732-4622	732-4624		
Purification of Serum Proteins, Enzymes									
	Affi-Gel Blue	Affinity	Cibacron Blue F3GA	0.2 ml serum 11 mg albumin	—	732-4642	732-4644		
Cartridge by Application	Support	Type of Separation	Chemical Form	Protein Capacity/ml*	Catalog Numbers by Cartridge Size				
					1 x 5 ml	5 x 5 ml	5 x 10 ml	1 x 50 ml	5 x 50 ml
Desalting, Buffer Exchange									
	Bio-Gel® P-6	Gel filtration	Polyacrylamide	10–22% CV**	732-4502	732-4504	732-5304	732-5312	732-5314

* Purification capacity based on bulk support; check individual instruction manuals for run conditions and specifications.

Specifications

Sizes (bed volumes)	1, 5, 10, and 50 ml
Maximum operating pressure	1, 5, and 10 ml: 45 psi 50 ml: 40 psi
Recommended flow rate	Varies for each support type*
Fittings	Female luer inlet and male luer outlet
Column material	Polypropylene
Frit material	Polyethylene (HDPE)
Dimensions (L x ID)**	1 ml — 40 mm x 5.6 mm 5 ml — 40 mm x 12.6 mm 10 ml — 80 mm x 12.6 mm*** 50 ml — 80 mm x 28.2 mm***

* Recommended flow rates are given for each support type in the instruction manual.

** L, length; ID, inner diameter.

*** Available for desalting cartridges only.

For current information on prepacked cartridges, visit www.bio-rad.com/cartridges/.

Cibacron is a trademark of Ciba-Geigy Corporation.

Purification and preparation of fusion proteins and affinity peptides containing at least two adjacent histidine residues may require a license under U.S. patent 5,284,933 and U.S. patent 5,310,663, including foreign patents (assignee: Hoffmann-La Roche).

Expression and purification of GST fusion proteins may require a license under U.S. patent 5,654,176 (assignee: Chemicon International).



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600 Canada 905 364 3435 China 86 20 8732 2339 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 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