

Get Your Fill.

A full line of empty chromatography columns and accessories.



Bio-Rad's Empty Column Line Is Packed With Options

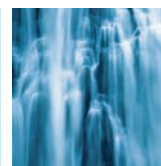
From low-pressure chromatography to high-resolution separations, we offer a comprehensive line of empty chromatography columns, accessories, and media to meet your specific chromatographic needs. As with all products from Bio-Rad, each is manufactured under the strictest quality control standards. You can expect only the highest quality with every purchase you make.

Empty Column Selection Guide

Application Requirement	Product Line	Pressure Range, psi	Volume Range, ml	Best Used With
High-resolution column	 Bio-Scale™ columns	0–1,000	1.9–22	High-performance system
General-purpose glass column	 Econo-Column® columns	0–14.7	0–1,374	Low-pressure or gravity-flow system
General-purpose disposable plastic column	 Econo-Pac® columns	0–14.7	0–20	Low-pressure or gravity-flow system
General-purpose disposable plastic column	 Poly-Prep® columns	0–14.7	0–2	Low-pressure or gravity-flow system
General-purpose disposable plastic column	 Bio-Spin® columns	—	0–1.2	Swinging bucket centrifuge
General-purpose disposable plastic column	 Micro Bio-Spin™ columns	—	0–0.8	Standard microcentrifuge

Bio-Rad also offers process-scale chromatography columns. Contact your local Bio-Rad office for more information.

Low-Pressure Chromatography Columns



Bio-Rad's Econo-Column line offers a complete range of versatile columns for low-pressure chromatography applications. A wide range of glass columns, jacketed columns, flow adaptors, reservoirs, and other chromatography essentials offers an unbeatable combination of performance, reliability, and value.

Glass Econo-Column Columns

Econo-Column chromatography columns are the standard in high-quality, affordable low-pressure chromatography columns.

- Available in lengths from 5 to 170 cm
- Inner diameter from 0.5 to 5.0 cm
- Volumes from 0 to 1,374 ml
- Operating pressure to 1 bar/14.7 psi
- Fully autoclavable and NaOH compatible

Jacketed Econo-Column Columns

Jacketed Econo-Column columns are ideal for chromatographic applications requiring temperature control, such as thermal chromatography of DNA separation using Bio-Gel® hydroxyapatite media.

- Integral water jacket for cooling or heating
- Open-ended column delivered with 2 flow adaptors
- Standard column delivered with 1 flow adaptor

Flow Adaptors

Flow adaptors significantly improve column performance by eliminating head space above the gel bed and by protecting the gel bed from disruption during sample loading.

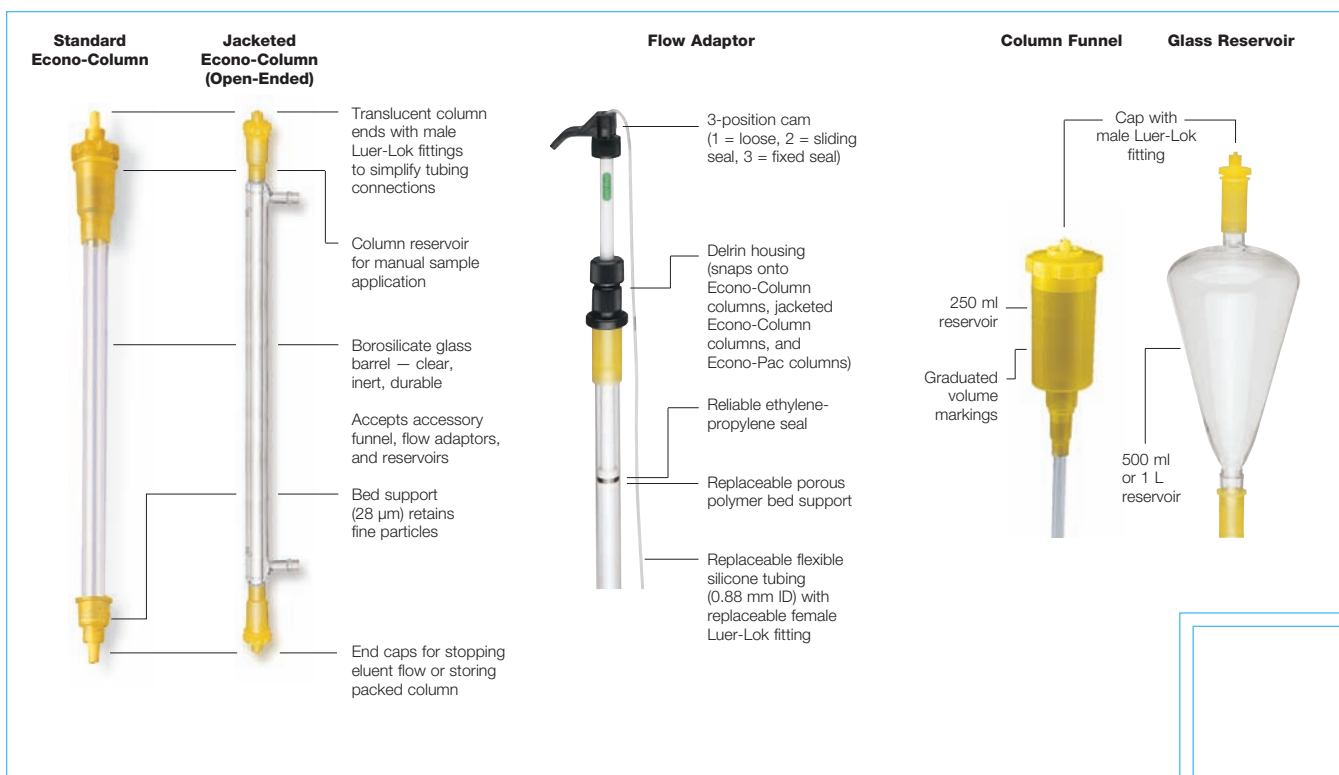
- Available for 1.0, 1.5, 2.5, and 5.0 cm columns
- Fit Econo-Column chromatography columns, jacketed Econo-Column chromatography columns, and Econo-Pac columns*

Column Funnel and Glass Reservoirs

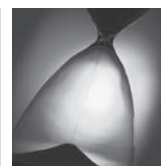
Econo-Column funnel and reservoirs are ideal for column packing, loading diluted samples, or delivering large volumes of buffer.

- Reservoirs available for 500 ml and 1 L capacities
- Reservoirs fit 0.5, 0.7, 1.0, and 1.5 cm ID Econo-Column columns
- Econo-Column funnel is ideal for packing columns up to 2.5 cm ID, including Econo-Column columns, Poly-Prep columns, and Econo-Pac columns

* Special Econo-Pac flow adaptor allows use with Econo-Pac columns.



Disposable Low-Pressure Chromatography Columns



All of Bio-Rad's disposable chromatography columns are made of chemically compatible polypropylene, have a 30 μm hydrophilic bed support, are fully autoclavable, and can be washed with NaOH. All are also available prepacked with selected media.

Bio-Spin and Micro Bio-Spin Columns

Bio-Spin and Micro Bio-Spin columns are disposable polypropylene spin columns that can be packed with a variety of chromatography media, such as gel filtration media, to remove dye terminators, clean up DNA primers, remove radiolabels, exchange buffers, or desalt proteins or DNA.

- Bio-Spin columns hold up to 1.2 ml of medium and fit in standard swinging bucket centrifuges
- Micro Bio-Spin columns hold up to 0.8 ml of medium and fit in standard microcentrifuges
- Both columns fit into standard microtubes
- Luer end fitting with snap-off tip
- Porous 30 μm polyethylene bed support retains fine particles
- Autoclavable and NaOH compatible

Poly-Prep Columns

Poly-Prep columns are convenient gravity-flow chromatography columns for sample preparation and small-scale chromatography applications, including work with radioisotopes and other applications where disposable products are required.

- 2 ml bed volume and integral 10 ml reservoir
- Porous 30 μm polyethylene bed support retains fine particles
- Luer end fitting with snap-off tip
- Autoclavable and NaOH compatible

Econo-Pac Columns

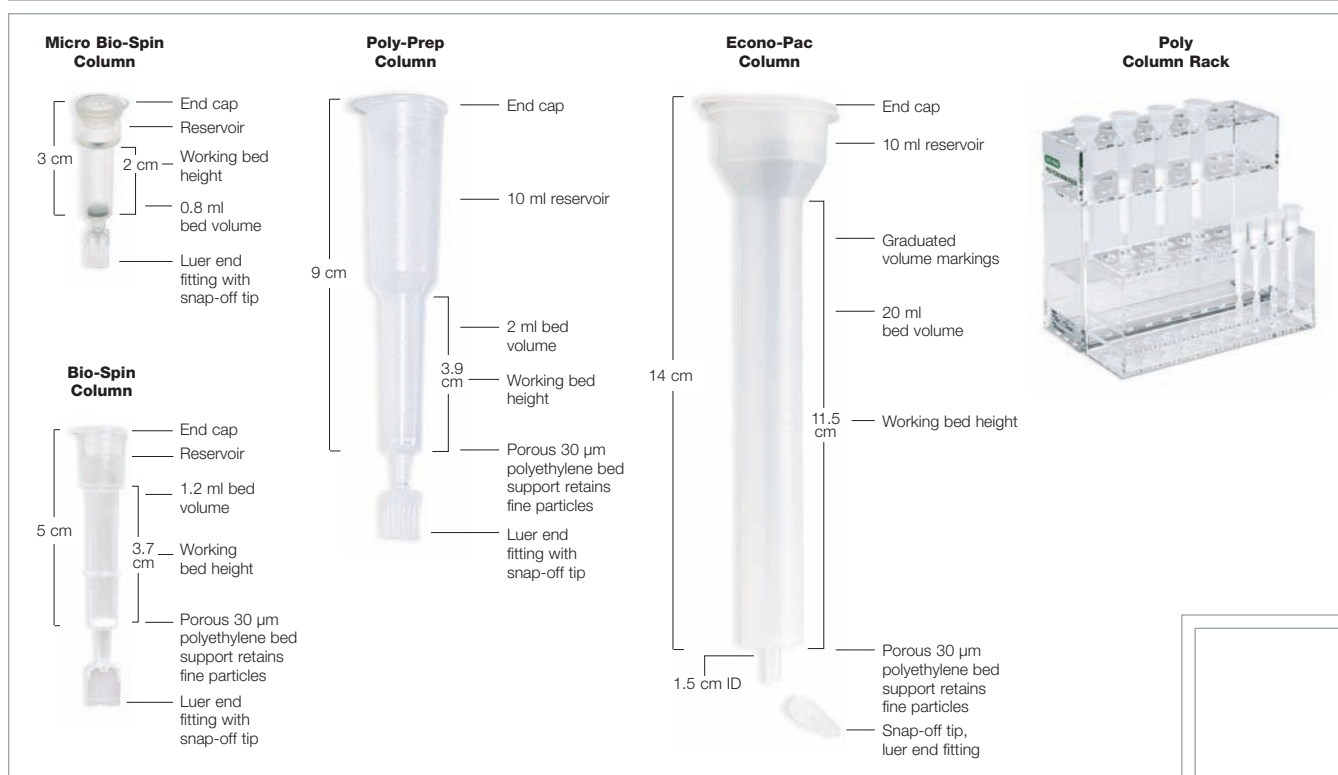
Econo-Pac columns are disposable 20 ml polypropylene columns for sample preparation and small-scale chromatography applications.

- 0–20 ml bed volume and 10 ml reservoir
- Porous 30 μm polyethylene bed support retains fine particles
- May be fitted with a flow adaptor or used for gravity-flow chromatography
- Luer end fitting with snap-off tip
- Autoclavable and NaOH compatible

Poly Column Rack

The poly column rack is ideal for using or storing Poly-Prep or Econo-Pac columns.

- Holds up to 20 Poly-Prep columns or 10 Econo-Pac columns
- Removable tube rack for Poly-Prep column use
- Reservoir tray to catch flow-through buffers



Medium-Pressure Chromatography Columns

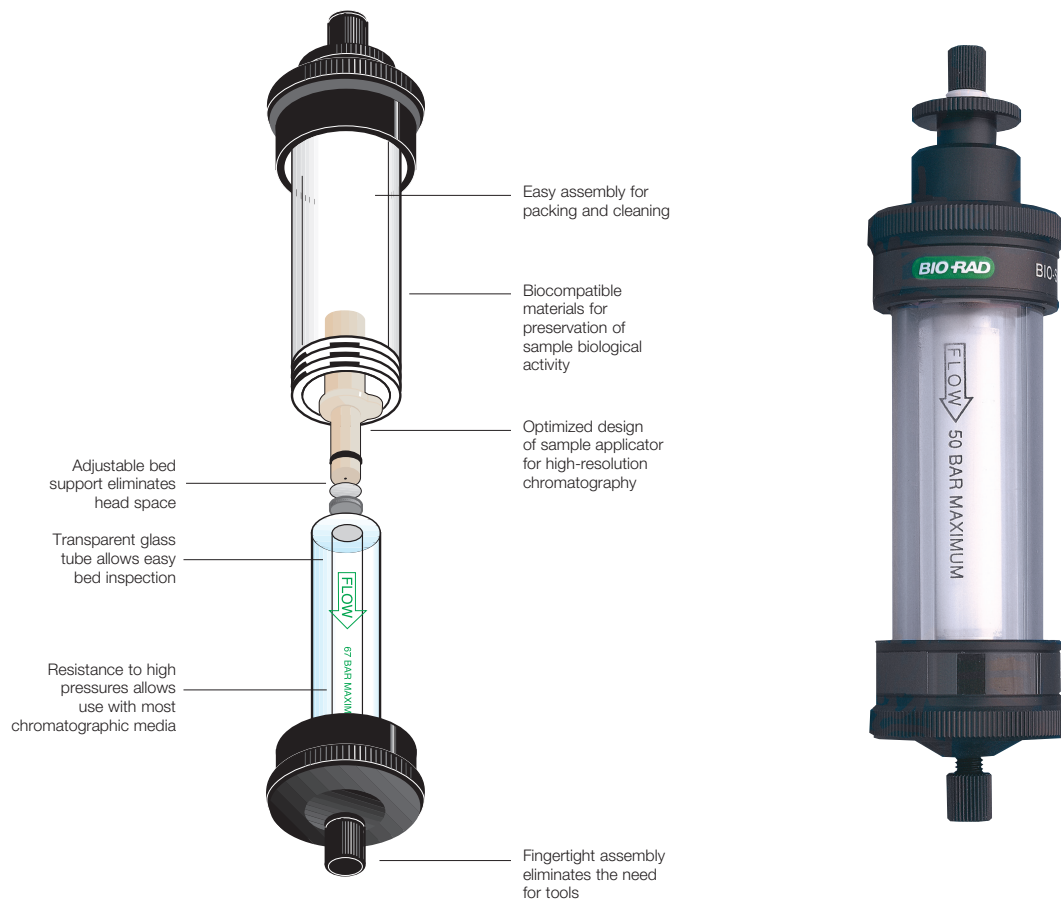


These high-quality columns are compatible with Bio-Rad's BioLogic™ systems or any medium- to high-pressure chromatography system. The range of columns has been specifically designed to provide high-resolution separation of biomolecules when packed with suitable chromatography media.

Bio-Scale MT High-Resolution Columns

Bio-Scale MT columns allow precise sample application and provide the low dead volume required for high-resolution separations.

- 4 column sizes (2, 5, 10, and 20 ml) allow easy scale-up of purification protocols
- Adjustable bed support eliminates head space
- Fingertight assembly eliminates the need for tools
- Biocompatible material resistant to high pressures allows use of most chromatographic media



Ordering Information

Catalog #	ID, cm	Length, cm	Cross-Sectional Area, cm ²	Maximum Volume, ml	Package Size
Econo-Column Chromatography Columns					
737-0507	0.5	5	0.20	1	2
737-0512	0.5	10	0.20	2	2
737-0517	0.5	15	0.20	3	2
737-0522	0.5	20	0.20	4	2
737-0707	0.7	5	0.39	2	2
737-0712	0.7	10	0.39	4	2
737-0717	0.7	15	0.39	6	2
737-0722	0.7	20	0.39	8	2
737-0732	0.7	30	0.39	12	2
737-0752	0.7	50	0.39	20	2
737-4506	0.5	5	0.20	1	4
737-4511	0.5	10	0.20	2	4
737-4516	0.5	15	0.20	3	4
737-4521	0.5	20	0.20	4	4
737-4706	0.7	5	0.39	2	4
737-4711	0.7	10	0.39	4	4
737-4716	0.7	15	0.39	6	4
737-4721	0.7	20	0.39	8	4
737-4731	0.7	30	0.39	12	4
737-4751	0.7	50	0.39	20	4
737-4006	1.0	5	0.79	4	4
737-1007	1.0	5	0.79	4	2
737-4011	1.0	10	0.79	8	4
737-1012	1.0	10	0.79	8	2
737-4021	1.0	20	0.79	16	4
737-1022	1.0	20	0.79	16	2
737-4031	1.0	30	0.79	24	4
737-1032	1.0	30	0.79	24	2
737-4051	1.0	50	0.79	40	4
737-1052	1.0	50	0.79	40	2
737-1091	1.0	100	0.79	79	2
737-1093	1.0	120	0.79	103	2
737-4150	1.5	5	1.77	9	4
737-1507	1.5	5	1.77	9	2
737-4151	1.5	10	1.77	18	4
737-1512	1.5	10	1.77	18	2
737-4156	1.5	15	1.77	27	4
737-1517	1.5	15	1.77	27	2
737-4152	1.5	20	1.77	35	4
737-1522	1.5	20	1.77	35	2
737-4153	1.5	30	1.77	53	4
737-1532	1.5	30	1.77	53	2
737-4155	1.5	50	1.77	89	4
737-1552	1.5	50	1.77	89	2
737-1576	1.5	75	1.77	124	2
737-1591	1.5	100	1.77	177	2
737-1593	1.5	120	1.77	230	2
737-1598	1.5	170	1.77	301	2
737-4250	2.5	5	4.91	25	4
737-2507	2.5	5	4.91	25	2
737-4251	2.5	10	4.91	49	4
737-2512	2.5	10	4.91	49	2
737-4252	2.5	20	4.91	98	4
737-2522	2.5	20	4.91	98	2
737-4253	2.5	30	4.91	147	4
737-2532	2.5	30	4.91	147	2
737-2551	2.5	50	4.91	246	2
737-2576	2.5	75	4.91	344	2
737-2591	2.5	100	4.91	491	2
737-2593	2.5	120	4.91	589	2
737-5011	5.0	10	19.63	196	1
737-5021	5.0	20	19.63	393	1
737-5031	5.0	30	19.63	589	1
737-5051	5.0	50	19.63	982	1
737-5071	5.0	70	19.63	1,374	1

Delrin is a trademark of E.I. du Pont de Nemours and Company. FPLC is a trademark of GE Healthcare. Luer-Lok is a trademark of Becton, Dickinson and Co.

Catalog #	Description	Volume, ml	Column Pressure Limit, psi		
Econo-Column Selection Packs					
737-6601	Econo-Column Selection Pack A, 1 each standard column, 0.7 x 10, 20, 30 cm; 1.5 x 30, 50 cm; 2.5 x 20, 50 cm				
737-6607	Econo-Column Selection Pack B, 1 each standard column, 1.0 x 20, 30, 50 cm; 1.5 x 20, 30, 50 cm				
Standard Econo-Column Jacketed Columns					
Catalog #	ID, cm	Length, cm	Cross-Sectional Area, cm ²	Maximum Volume, ml	Package Size
737-6108	0.7	15	0.37	6	1
737-6116	1.0	15	0.79	12	1
737-6131	1.0	30	0.79	25	1
737-6151	1.5	50	1.77	89	1
Econo-Column Open-Ended Jacketed Column*					
737-6201	1.0	30	0.79	25	1
Catalog #	Description	Column ID, cm	Functional Length, cm		
Flow Adaptors					
738-0014	Flow Adaptor	1.0	1-7		
738-0015	Flow Adaptor	1.0	1-14		
738-0016	Flow Adaptor	1.5	1-14		
738-0017	Flow Adaptor	2.5	1-14		
738-0018	Flow Adaptor**	5.0	1-14		
738-0019	Econo-Pac Flow Adaptor	1.5	—		
Flow Adaptor Replacement Parts					
738-0022	Flow Adaptor Maintenance Kit, for 5.0 cm flow adaptor, includes 2 bed supports, 2 O-rings				
738-0024	Flow Adaptor Maintenance Kit, for 1.0 cm adaptor with cam mechanism, includes 10 bed supports, 5 O-rings				
738-0025	Flow Adaptor Maintenance Kit, for 1.5 cm adaptor with cam mechanism, includes 10 bed supports, 5 O-rings				
738-0027	Flow Adaptor Maintenance Kit, for 2.5 cm adaptor with cam mechanism, includes 10 bed supports, 5 O-rings				
Funnel and Reservoirs					
731-0003	Econo-Column Funnels, 250 ml, 5				
737-9112	Econo-Column Reservoir, 500 ml				
737-9113	Econo-Column Reservoir, 1 L				
Catalog #	Description	Volume, ml	Column Pressure Limit, psi		
Bio-Scale MT Columns					
751-0081	Bio-Scale MT2 Column, 7 x 52 mm	1.9-2.3	1,000		
751-0083	Bio-Scale MT5 Column, 10 x 64 mm	4.6-5.7	750		
751-0085	Bio-Scale MT10 Column, 12 x 88 mm	9.5-11.3	600		
751-0087	Bio-Scale MT20 Column, 15 x 113 mm	19.4-21.9	500		
Bio-Scale Kits					
751-0091	Bio-Scale 2 Replacement Parts Kit, includes 5 frits, 5 distribution screens, 2 O-rings, 1 frit remover				
751-0093	Bio-Scale 5 Replacement Parts Kit				
751-0095	Bio-Scale 10 Replacement Parts Kit				
751-0097	Bio-Scale 20 Replacement Parts Kit				
750-0554	1/16" OD (1.6 mm) Post-Pump Fittings, includes Delrin nut, ferrules, lock ring, 10 sets				
750-0568	UNO® 10-32 Fittings Kit, set of 2 nuts and 4 ferrules to connect a Bio-Scale MT column to an HPLC system				
750-0567	UNO M6 Fittings Kit, set of 2 nuts and 4 ferrules to connect a Bio-Scale MT column to an FPLC system				
Bio-Spin and Micro Bio-Spin Columns					
732-6008	Bio-Spin Chromatography Columns, empty, 100				
732-6025	Bio-Spin Chromatography Columns, empty, 1,000				
732-6204	Micro Bio-Spin Chromatography Columns, empty, 100				
731-1660	End Caps, for Micro Bio-Spin chromatography columns, 1,000				
Econo-Pac Columns					
732-1010	Econo-Pac Chromatography Columns, empty, includes upper bed supports, end caps, tip closures, 50				
732-1011	Econo-Pac Chromatography Columns, empty, includes upper bed supports, end caps, tip closures, 500				
Poly-Prep Columns					
731-1550	Poly-Prep Chromatography Columns, empty, includes end caps and tip closures, 50				
731-1553	Poly-Prep Chromatography Columns, empty, includes end caps and tip closures, 1,000				
Accessories					
731-1555	Poly-Prep Column Stack Cap, 50				
731-7005	Poly Column Rack, 20 place, with removable tube rack				
732-8102	2-Way Stopcock, 10				
731-8232	Female Luer Plug, 25				

* Includes 2 flow adaptors.

** 5.0 cm flow adaptor does not include cam mechanism.



**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