

# Econo Alpha Empty Chromatography Columns

**Pack your choice of chromatography resin in a column size that fits your needs.**

Bio-Rad offers a wide assortment of resins and prepacked low- and medium-pressure columns for affinity, ion exchange, size exclusion, mixed-mode, and hydrophobic interaction chromatography. With Econo Alpha Chromatography Columns you can customize your low- to medium-pressure chromatography applications.

## Features and Benefits

- Robust and easy-to-pack hardware
- Compatible with common aqueous buffers used in protein purification
- Wide range of accessories available for order

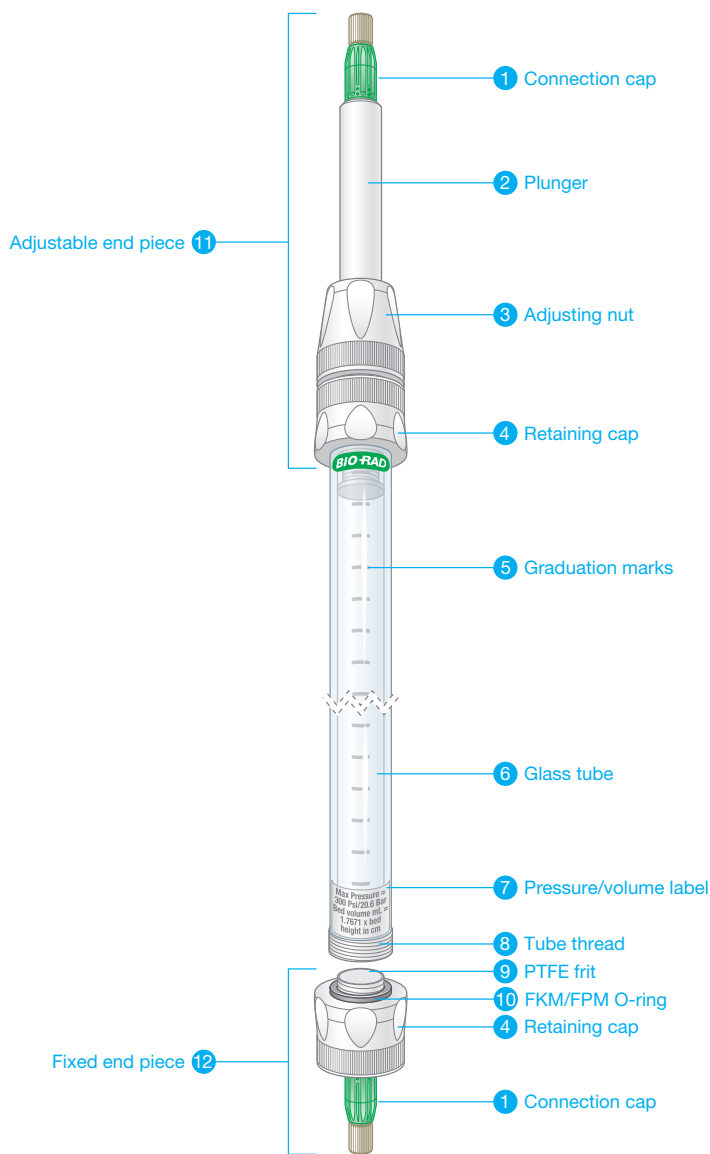
## Applications and Uses

- Low- to medium-pressure chromatography
- Designed to work seamlessly with the NGC Chromatography System family
- Columns come with preinstalled 10–32 and 1/4–28 fittings for maximum compatibility with other fast protein liquid chromatography systems

## Standard Glass Column Configuration

Econo Alpha Chromatography Columns have one fixed end piece and one adjustable end piece (plunger), to adjust bed height, and frits with a 5 µm pore size.





Econo Alpha Column

**Specifications**

Parameter	Description
Operating temperature	4–40°C
pH stability	1–14
Chemical stability	Resistant to aqueous solutions, including 8 M urea and 6 M guanidine hydrochloride, and most solvents used in liquid chromatography. Not resistant to acetone, ketones, chlorinated hydrocarbons, aliphatic ester, phenol, >10% NaOH, >10% HCl, >5% acetic acid, or strong mineral acid

**Ordering Information**

**Assembled Columns**

Catalog Number	Column, mm	Bed Height, cm		Bed Volume, ml		Pressure Rating	
		Min	Max	Min	Max	psi	MPa
12009463	6.6 x 50	0.24	2	0.1	0.7	900	6.20
12009429	10 x 100	0.2	7	0.2	5.5	600	4.13
12009430	15 x 100	0.2	7	0.4	12.4	300	2.06
12009461	25 x 100	0.2	7	1.2	34.4	150	1.03
12009462	10 x 500	38.8	47	30.6	36.9	600	4.13
12009427	15 x 500	38.8	47	68.9	83.1	300	2.06
12009428	25 x 500	38.8	47	191.4	230.7	150	1.03

**Accessories**

Column packing accessories, additional plungers, frits (5–50 µm pore size), and O-rings.



Visit [bio-rad.com/Lab-Scale](http://bio-rad.com/Lab-Scale) or [bio-rad.com/Process-Scale](http://bio-rad.com/Process-Scale) to view our resin offerings.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

All trademarks used herein are the property of their respective owner.



**Bio-Rad Laboratories, Inc.**

Life Science Group

Website [bio-rad.com](http://bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

