

Econo Alpha Empty Glass Chromatography Columns

Catalog #	Description
Multiple	Econo Alpha Empty Glass Chromatography Columns, multiple sizes

For research purposes only.

Overview

Econo Alpha Columns are convenient, easy-to-pack chromatography columns. They are supplied assembled and ready for use. Refer to the instruction manual (10000108068) and go to bio-rad.com/ResinsandColumns for additional information.

Chemical Stability

Econo Alpha Columns are resistant to aqueous solutions, including 8 M urea and 6 M guanidine hydrochloride, and most solvents used in liquid chromatography. These columns are not resistant to acetone, ketones, chlorinated hydrocarbons, aliphatic ester, phenol, >10% NaOH, >10% HCl, >5% acetic acid, or strong mineral acids.

Specifications

Property	Operating Parameter
Operating temperature	4–40°C
pH stability	1–14
Component	Composition
Glass column	Borosilicate glass
End piece and plunger	Polytetrafluoroethylene (PTFE)
Frit (bed support; 5 µm)	PTFE
O-ring	Fluorocarbon/fluoropolymer
Adjusting and retaining nuts	Acetal
Column Size	Operating Pressure
6.6 mm	900 psi, 62.05 bar, 6.20 MPa
10 mm	600 psi, 41.36 bar, 4.13 MPa
15 mm	300 psi, 20.68 bar, 2.06 MPa
25 mm	150 psi, 10.34 bar, 1.03 MPa

Configuration

The Econo Alpha Column is supplied assembled and ready for use, with fittings to connect the column into an existing system. It ships in a configuration that consists of one fixed end piece as a bed support and one adjustable end piece (plunger) to adjust the bed height. The adjustable plunger allows precise adjustment with minimal disturbance to the packed bed. Additional adjustable end pieces may be purchased separately to enable even more flexibility in bed height.

Column Bed Heights and Volumes

Catalog Number	Column, mm	Bed Height, cm		Bed Volume, ml	
		Minimum	Maximum	Minimum	Maximum
12009463	6.6 x 50	0.24	2	0.1	0.7
12009429	10 x 100	0.2	7	0.2	5.5
12009430	15 x 100	0.2	7	0.4	12.4
12009461	25 x 100	0.2	7	1.2	34.4
12009462	10 x 500	38.8	47	30.6	36.9
12009427	15 x 500	38.8	47	68.9	83.1
12009428	25 x 500	38.8	47	191.4	230.7

Ordering Information

Catalog #	Description
Econo Alpha Columns	
12009463	Econo Alpha Column, 6.6 x 50 mm
12009429	Econo Alpha Column, 10 x 100 mm
12009430	Econo Alpha Column, 15 x 100 mm
12009461	Econo Alpha Column, 25 x 100 mm
12009462	Econo Alpha Column, 10 x 500 mm
12009427	Econo Alpha Column, 15 x 500 mm
12009428	Econo Alpha Column, 25 x 500 mm
Adjustable Plungers	
12009493	Econo Alpha Adjustable Plunger, 10 mm
12009492	Econo Alpha Adjustable Plunger, 15 mm
12009494	Econo Alpha Adjustable Plunger, 25 mm
Packing Sleeves (connector)	
12009487	Econo Alpha Packing Sleeve, 6.6 mm
12009486	Econo Alpha Packing Sleeve, 10 mm
12009484	Econo Alpha Packing Sleeve, 15 mm
12009485	Econo Alpha Packing Sleeve, 25 mm
Column Extensions	
12009501	Econo Alpha Column Extension, 6.6 x 50 mm
12009497	Econo Alpha Column Extension, 10 x 100 mm
12009498	Econo Alpha Column Extension, 15 x 100 mm
12009499	Econo Alpha Column Extension, 25 x 100 mm
12009500	Econo Alpha Column Extension, 10 x 500 mm
12009495	Econo Alpha Column Extension, 15 x 500 mm
12009496	Econo Alpha Column Extension, 25 x 500 mm
Frits	
12009479	Econo Alpha Frit, 6.6 mm, 5 µm
12009480	Econo Alpha Frit, 6.6 mm, 10 µm
12009481	Econo Alpha Frit, 6.6 mm, 20 µm
12009482	Econo Alpha Frit, 6.6 mm, 30 µm
12009483	Econo Alpha Frit, 6.6 mm, 50 µm
12009474	Econo Alpha Frit, 10 mm, 5 µm
12009475	Econo Alpha Frit, 10 mm, 10 µm
12009476	Econo Alpha Frit, 10 mm, 20 µm
12009477	Econo Alpha Frit, 10 mm, 30 µm
12009478	Econo Alpha Frit, 10 mm, 50 µm
12009464	Econo Alpha Frit, 15 mm, 5 µm
12009465	Econo Alpha Frit, 15 mm, 10 µm
12009466	Econo Alpha Frit, 15 mm, 20 µm
12009467	Econo Alpha Frit, 15 mm, 30 µm
12009468	Econo Alpha Frit, 15 mm, 50 µm
12009469	Econo Alpha Frit, 25 mm, 5 µm
12009470	Econo Alpha Frit, 25 mm, 10 µm
12009471	Econo Alpha Frit, 25 mm, 20 µm
12009472	Econo Alpha Frit, 25 mm, 30 µm
12009473	Econo Alpha Frit, 25 mm, 50 µm
O-Rings	
12009491	Econo Alpha O-Rings, 6.6 mm
12009490	Econo Alpha O-Rings, 10 mm
12009488	Econo Alpha O-Rings, 15 mm
12009489	Econo Alpha O-Rings, 25 mm

Visit bio-rad.com/ResinsandColumns for the most up-to-date list of column and resin offerings.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.
All trademarks used herein are the property of their respective owner.