

Bio-Rad, Affi-Gel, Affi-Prep, Bio-Gel, CHT, and Macro-Prep are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. Nuvia aPrime 4A Resin is covered by U.S. Patent Number 9,669,402 and foreign counterparts. All trademarks used herein are the property of their respective owner.

Resin	Chemistry	Media	Chemical Form	Protein Capacity	EconoFit Catalog Numbers by Column Size	
Mixed mode	Nuvia aPrime 4A	40 µm	Hydrophobic anion	At 300 cm/hr: >50 mg acidic mAb/ml	12009280	
				At 300 cm/hr: >40 mg human IgG/ml	12009281	
	MPC I, 40 µm	40 µm	Ca ²⁺ , PO ₄ ⁺ , OH, F	≥25 mg lysozyme/g; 25–60 mg IgG/g	12009279	
				14–21.5 mg/g	12009252	
	CFT Type II, 80 µm	40 µm	Ca ₁₀ (PO ₄) ₆ F ₂	Lysozyme: 15–25 mg IgG/g	12009260	
				≥12.5 mg lysozyme/g; 25–60 mg IgG/g	12009256	
	CFT Type I, 80 µm	40 µm	Ca ²⁺ , PO ₄ ⁺ , OH	≥25 mg lysozyme/g; 17–25 mg lysozyme/g	12009261	
				At 500 cm/hr: mAb G/ml (pI 9.2) At 100 cm/hr: >60 mg human IgG/ml	12009259	
	CFT Type II, 40 µm	40 µm	Ca ²⁺ , PO ₄ ⁺ , OH	15–25 mg human IgG/ml	12009257	
				25–60 mg human IgG/ml	12009254	
					1 x 1 ml	5 x 1 ml
					1 x 5 ml	5 x 5 ml

Ordering Information



EconoFit Chromatography Columns

Catalog #	Description
Multiple	EconoFit Chromatography Columns, 1 and 5 ml

For research purposes only.

Overview

EconoFit Chromatography Columns are convenient, disposable, prepacked low-pressure chromatography columns that offer increased run-to-run uniformity through the column design. This product insert will guide you in the basic use of the columns. For additional information, including resin-specific guidance and the most up-to-date list of bottled resin and column offerings, go to bio-rad.com/ResinsandColumns and consult the instruction manual. Larger package sizes of media are available for process-scale chromatography.

Specifications

Property	Description
Dimensions (D x H)	1 ml: 7 x 25 mm 5 ml: 16 x 25 mm
Maximum pressure tolerance (hardware)	72 psi/5 bar/0.5 MPa
Recommended flow rate	Varies for each support type*
Fittings	10-32 (1/16"), female inlet and male outlet
Column material	Polypropylene
Frit material	High-density polyethylene

* Please refer to the online product manual for resin-specific guidance.

Ordering Information

Resin Chemistry	Media	Chemical Form	Protein Capacity	EconoFit Catalog Numbers by Column Size			
				1 x 1 ml	5 x 1 ml	1 x 5 ml	5 x 5 ml
Affinity	Affi-Gel Blue	Cibacron Blue F3GA	0.2 ml serum, 11 mg albumin	—	—	12009234	12009235
	Affi-Prep Protein A	Protein A	8–10 mg mouse monoclonal IgG1	—	12009236	12009237	—
	DEAE Affi-Gel Blue	DEAE–Cibacron Blue	0.2–1.0 ml serum F3GA and DEAE	—	—	12009262	12009263
	Nuvia IMAC Ni-Charged	Nitrilotriacetic acid, Ni-charged	≥40 mg/ml recombinant histidine-tagged protein	12009288	12009285	12009286	12009287
	Nuvia IMAC Uncharged	Nitrilotriacetic acid	N/A	12009289	—	—	—
	Profinity eXact	Modified subtilisin protease	>3 mg/ml tag-free maltose binding protein	12009292	12009293	12009294	—
	Profinity GST	Glutathione	≥7 mg/ml GST-tagged protein	—	12009295	12009296	12009297
	Profinity IMAC	Iminodiacetic acid	≥15 mg/ml recombinant histidine-tagged protein	—	12009298	12009299	12009300
	UNOsphere SUPrA	rProtein A	25–30 mg IgG/ml	—	12009322	12009323	—
Ion exchange	Macro-Prep DEAE	–NH ⁺ (C ₂ H ₅) ₂	>35 mg BSA	12009274	12009264	12009265	12009266
	Macro-Prep High Q	–N ⁺ (CH ₃) ₃	≥37 mg BSA/ml	12009275	12009267	12009268	12009269
	Macro-Prep High S	–SO ₃ [–]	≥49 mg human IgG	12009276	12009270	12009271	12009272
	Macro-Prep High Q-3HT	–N ⁺ (CH ₃) ₃	≥37 mg BSA/ml	12009283	—	—	—
	UNOsphere Q	–N ⁺ (CH ₃) ₃	At 150 cm/hr: 180 mg BSA/ml At 300 cm/hr: 125 mg BSA/ml	12009307	12009301	12009302	12009303
	UNOsphere S	–SO ₃ [–]	At 150 cm/hr: 60 mg IgG/ml At 300 cm/hr: 30 mg IgG/ml	12009308	12009304	12009305	12009306
	Nuvia Q	–N ⁺ (CH ₃) ₃	At 300 cm/hr: >170 mg BSA/ml	12009290	—	—	—
	Nuvia S	–SO ₃ [–]	At 300 cm/hr: >110 mg human IgG/ml	12009291	—	—	—
	Nuvia HR-S	–SO ₃ [–]	At 300 cm/hr: ≥70 mg human IgG/ml	12009284	—	—	—
	Nuvia HP-Q	–N ⁺ (CH ₃) ₃	At 100 cm/hr: >50 mg thyroglobulin/ml	12009282	—	—	—
Hydrophobic interaction	Macro-Prep Methyl	–CH ₃	>25 mg HSA/ml	12009277	—	—	—
	Macro-Prep t-Butyl	–C(CH ₃) ₃	>15 mg HSA/ml	12009278	—	—	—
Desalting/size exclusion	Bio-Gel P-6 Desalting	Polyacrylamide	10–22% column volume	—	—	12009238	12009239
Empty				—	12009232	—	12009233

continues