



Resin and Column Selection Guide

The right column for the right resin



Resin and Column Selection Guide

Affinity | Desalting/Buffer Exchange



	EconoFit Columns	Foresight Columns*	Econo Alpha Columns	Bio-Scale MT Columns
Main features	Prepacked and empty columns for automated or manual chromatography	Robust prepacked columns for high-throughput experiments	Low- to medium-pressure chromatography; glass; adjustable bed height	High-resolution chromatography; acrylic jacketed glass; adjustable bed height
Bed volume	1/5 ml	1/5 ml	0.1–230.7 ml	1.9–21.9 ml
Bed height	2.5 cm	2/10 cm	0.2–47 cm	5–12.4 cm
Maximum operating pressure	72 psi/0.5 MPa	290–435 psi/2–3 MPa	150–900 psi/ 1.03–6.20 MPa	500–1,000 psi/ 3.45–6.9 MPa
Frit pore size	15 µm	7–12 µm	5–50 µm	10 µm
Adaptors available for column	Yes	Yes	Yes	Yes
Available prepacked	Yes	Yes	No	Yes

Chromatography Technique

Affinity Stationary Phase

Nuvia IMAC	P	P	R	N
Nuvia IMAC, Uncharged	P	—	R	N
Profinity IMAC	P	—	R	N
Profinity IMAC, Uncharged	R	—	R	N
Profinity GST**	P	—	N	N
Profinity eXact	P	—	R	N
Profinity Epoxide	R	—	R	R
UNOsphere SUPrA	P	—	R	N
Affi-Gel Protein A	N	—	N	N
Affi-Prep Protein A	P	—	R	R
DEAE Affi-Gel Blue	P	—	R	N
CM Affi-Gel Blue	R	—	R	N
Affi-Gel Blue	P	—	R	N
Affi-Gel Boronate	R	—	R	N
Affi-Prep Polymyxin	R	—	R	N
Affi-Gel Heparin	R	—	R	N
Affi-Gel Hz Hydrazide	R	—	R	N
Affi-Gel 102	R	—	R	N
Affi-Gel 10/15	R	—	R	N

Desalting/Buffer Exchange Stationary Phase

Bio-Gel P-6 Desalting***	P	—	R	N
Bio-Gel P-30	R	—	R	N

* Foresight Columns are available only in prepacked format.

** Available only in prepacked format.

*** Also available in Bio-Scale Mini 10 and 50 ml prepacked formats.

P, available prepacked; **R**, recommended; **N**, not recommended; **—**, empty column not available

Resin and Column Selection Guide

Affinity | Desalting/Buffer Exchange



	Micro/Mini/Bio-Spin Columns	Poly-Prep and Econo-Pac Columns	Glass Econo-Column Columns
Main features	Broad range of sizes in an economical format for use in a standard centrifuge	Empty Econo-Pac and Poly-Prep Gravity Flow Columns are ideal for sample and buffer preparation and for gravity flow chromatography	Broad range of sizes in an economical format with a porous polymer bed support
Maximum bed volume	0.8/1.2/1.6 ml	2/20 ml	1,374 ml
Bed height	2/2.8/3.7 cm	3.9/11.5 cm	5–170 cm
Maximum operating pressure	1,000 x g spin	14.7 psi/0.1 MPa	14.7 psi/0.1 MPa
Frit pore size	30 µm	30 µm	30 µm
Adapters available for column	N/A	Yes	Yes
Available prepacked	Yes	Yes	No

Chromatography Technique

Affinity Stationary Phase

Nuvia IMAC	R	R	R
Nuvia IMAC, Uncharged	R	R	R
Profinity IMAC	N	R	R
Profinity IMAC, Uncharged	N	R	R
Profinity GST*	N	N	N
Profinity eXact	P	R	R
Profinity Epoxide	N	R	R
UNOsphere SUPrA	R	R	R
Affi-Gel Protein A	R	P	R
Affi-Prep Protein A	N	N	N
DEAE Affi-Gel Blue	R	R	R
CM Affi-Gel Blue	R	R	R
Affi-Gel Blue	R	R	R
Affi-Gel Boronate	R	R	R
Affi-Prep Polymyxin	R	R	R
Affi-Gel Heparin	R	R	R
Affi-Gel Hz Hydrazide	R	R	R
Affi-Gel 102	R	R	R
Affi-Gel 10/15	R	R	R

Desalting/Buffer Exchange Stationary Phase

Bio-Gel P-6 Desalting	P	P	R
Bio-Gel P-30	P	R	R

* Available only in prepacked format.

P, available prepacked; R, recommended; N, not recommended

Resin and Column Selection Guide

Ion Exchange | Size Exclusion/Gel Filtration



	EconoFit Columns	Foresight Columns*	Econo Alpha Columns	Bio-Scale MT Columns
Main features	Prepacked and empty columns for automated or manual chromatography	Robust prepacked columns for high-throughput experiments	Low- to medium-pressure chromatography; glass; adjustable bed height	High-resolution chromatography; acrylic jacketed glass; adjustable bed height
Bed volume	1/5 ml	1/5 ml	0.1–230.7 ml	1.9–21.9 ml
Bed height	2.5 cm	2/10 cm	0.2–47 cm	5–12.4 cm
Maximum operating pressure	72 psi/0.5 MPa	290–435 psi/2–3 MPa	150–900 psi/1.03–6.20 MPa	500–1,000 psi/3.45–6.9 MPa
Frit pore size	15 µm	7–12 µm	5–50 µm	10 µm
Adaptors available for column	Yes	Yes	Yes	Yes
Available prepacked	Yes	Yes	No	Yes

Chromatography Technique

Ion Exchange Stationary Phase

Macro-Prep High Q	P	–	R	N
Macro-Prep High Q-3HT	P	–	R	N
Macro-Prep DEAE	P	–	R	N
Macro-Prep High S	P	–	R	N
Macro-Prep CM	P	–	R	N
Macro-Prep 25 Q	N	–	R	R
Macro-Prep 25 S	N	–	R	R
Nuvia Q	P	P	R	N
Nuvia HP-Q	P	P	R	N
Nuvia S	P	P	R	N
Nuvia HR-S	P	P	R	N
UNOsphere Q	P	P	R	N
UNOsphere S	P	P	R	N
AG 1	N	–	R	N
AG 2	N	–	R	N
AG 50W	N	–	R	N
AG MP-50	N	–	R	N
AG 501	N	–	R	N
Bio-Rex	N	–	R	N
Chelex	N	–	R	N
Bio-Rex 70	N	–	R	N
Bio-Rex RG 501-X8	N	–	R	N

Size Exclusion/Gel Filtration Stationary Phase

Bio-Gel P	P	–	R	N
Bio-Gel A	R	–	R	N
Bio-Beads S-X	R	–	R	N

Other Prepacked Columns*

ENrich Q and S	High-resolution ion exchange chromatography; bed volume: 1/8 ml; maximum operating pressure: 500 psi/3.45 MPa
ENrich SEC 70 and 650	High-resolution size exclusion chromatography; bed volume: 24 ml; maximum operating pressure: 600 psi/4.1 MPa
UNO Q and S	Continuous bed ion exchange chromatography; bed volume: 1.3/6/12 ml; maximum operating pressure: 700 psi/4.82 MPa
UNO Q and S Polishing	Continuous bed ion exchange chromatography; bed volume: 0.16 ml; maximum operating pressure: 200 psi/1.3 MPa

* Foresight, ENrich, and UNO Columns are available only in prepacked format.

P, available prepacked; **R**, recommended; **N**, not recommended; **–**, empty column not available

Resin and Column Selection Guide

Ion Exchange | Size Exclusion/Gel Filtration



	Micro/Mini/Bio-Spin Columns	Poly-Prep and Econo-Pac Columns	Glass Econo-Column Columns
Main features	Broad range of sizes in an economical format for use in a standard centrifuge	Empty Econo-Pac and Poly-Prep Gravity Flow Columns are ideal for sample and buffer preparation and for gravity flow chromatography	Broad range of sizes in an economical format with a porous polymer bed support
Bed volume	0.8/1.2/1.6 ml	2/20 ml	1,374 ml
Bed height	2/2.8/3.7 cm	3.9/11.5 cm	5–170 cm
Maximum operating pressure	1,000 x g spin	14.7 psi/0.1 MPa	14.7 psi/0.1 MPa
Frit pore size	30 µm	30 µm	30 µm
Adapters available for column	N/A	Yes	Yes
Available prepacked	Yes	Yes	No

Chromatography Technique

Ion Exchange Stationary Phase

Macro-Prep High Q	R	R	R
Macro-Prep High Q-3HT	R	R	R
Macro-Prep DEAE	R	R	R
Macro-Prep High S	R	R	R
Macro-Prep CM	R	R	R
Macro-Prep 25 Q	N	N	N
Macro-Prep 25 S	N	N	N
Nuvia Q	R	R	R
Nuvia HP-Q	R	R	R
Nuvia S	R	R	R
Nuvia HR-S	R	R	R
UNOsphere Q	R	R	R
UNOsphere S	R	R	R
AG 1	R	P	R
AG 2	R	R	R
AG 50W	R	P	R
AG MP-50	R	R	R
AG 501	R	R	R
Bio-Rex	R	R	R
Chelex	R	R	R
Bio-Rex 70	N	R	R
Bio-Rex RG 501-X8	N	R	R

Size Exclusion/Gel Filtration Stationary Phase

Bio-Gel P	P	P	R
Bio-Gel A	R	R	R
Bio-Beads S-X	R	R	R

P, available prepacked; R, recommended; N, not recommended

Resin and Column Selection Guide

Mixed-Mode | Hydrophobic Interaction



	EconoFit Columns	Foresight Columns*	Econo Alpha Columns	Bio-Scale MT Columns
Main features	Prepacked and empty columns for automated or manual chromatography	Robust prepacked columns, for high-throughput experiments	Low- to medium-pressure chromatography; glass; adjustable bed height	High-resolution chromatography; acrylic jacketed glass; adjustable bed height
Bed volume	1/5 ml	1/5 ml	0.1–230.7 ml	1.9–21.9 ml
Bed height	2.5 cm	2/10 cm	0.2–47 cm	5–12.4 cm
Maximum operating pressure	72 psi/0.5 MPa	290–435 psi/2–3 MPa	150–900 psi/ 1.03–6.20 MPa	500–1,000 psi/ 3.45–6.9 MPa
Frit pore size	15 µm	7–12 µm	5–50 µm	10 µm
Adapters available for column	Yes	Yes	Yes	Yes
Available prepacked	Yes	Yes	No	Yes**

Chromatography Technique

Mixed-Mode Stationary Phase

CHT Type I, 20 µm	N	–	R	R
CHT Type II, 20 µm	N	–	R	R
CHT Type I, 40 or 80 µm	P	P	R	R
CHT Type II, 40 or 80 µm	P	P	R	R
CHT XT, 40 µm	P	P	R	R
MPC Type I, 40 µm	P	P	R	R
CFT Type II, 40 µm	P	–	R	R
Nuvia aPrime 4A	P	P	R	N
Nuvia cPrime	P	P	R	N
Bio-Gel HT	N	–	R	N
Bio-Gel HTP	N	–	R	N

Hydrophobic Interaction Stationary Phase

Macro-Prep t-Butyl	P	–	R	N
Macro-Prep Methyl	P	–	R	N
Bio-Beads SM-2	R	–	R	N

* Foresight Columns are available only in prepacked format.

** Available prepacked only with CHT Type I, 10 µm stationary phase. See ordering information at the end of this document.

P, available prepacked; **R**, recommended; **N**, not recommended; **–**, empty column not available

Resin and Column Selection Guide

Mixed-Mode | Hydrophobic Interaction



	Micro/Mini/Bio-Spin Columns	Poly-Prep and Econo-Pac Columns	Glass Econo-Column Columns
Main features	Broad range of sizes in an economical format for use in a standard centrifuge	Empty Econo-Pac and Poly-Prep Gravity Flow Columns are ideal for sample and buffer preparation and for gravity flow chromatography	Broad range of sizes in an economical format with a porous polymer bed support
Bed volume	0.8/1.2/1.6 ml	2/20 ml	1,374 ml
Bed height	2/2.8/3.7 cm	3.9/11.5 cm	5–170 cm
Maximum operating pressure	1,000 x g spin	14.7 psi/0.1 MPa	14.7 psi/0.1 MPa
Frit pore size	30 µm	30 µm	30 µm
Adapters available for column	N/A	Yes	Yes
Available prepacked	Yes	Yes	No

Chromatography Technique

Mixed-Mode Stationary Phase

CHT Type I, 20 µm	N	N	N
CHT Type II, 20 µm	N	N	N
CHT Type I, 40 or 80 µm	N	R	R
CHT Type II, 40 or 80 µm	N	R	R
CHT XT, 40 µm	N	R	R
MPC Type I, 40 µm	N	R	R
CFT Type II, 40 µm	N	R	R
Nuvia aPrime 4A	N	N	R
Nuvia cPrime	N	N	R
Bio-Gel HT	N	R	R
Bio-Gel HTP	N	R	R

Hydrophobic Interaction Stationary Phase

Macro-Prep t-Butyl	R	R	R
Macro-Prep Methyl	R	R	R
Bio-Beads SM-2	R	R	R

P, available prepacked; R, recommended; N, not recommended

Ordering Information

Description	Catalog #	Quantity
Prepacked Columns		
Gel Filtration/Desalting/Small Molecule Removal		
ENrich SEC 70	7801070	1 x 24 ml
ENrich SEC 650	7801650	1 x 24 ml
EconoFit Bio-Gel P-6 Desalting	12009238 12009239	1 x 5 ml 5 x 5 ml
Bio-Scale Mini Bio-Gel P-6 Desalting	7325304 7325312 7325314	5 x 10 ml 1 x 50 ml 5 x 50 ml
Econo-Pac 10DG Desalting	7322010	30 x 10 ml
Micro Bio-Spin P-6 Gel, Tris	7326221 7326222 7326225	25 x 0.8 ml 100 x 0.8 ml 1,000 x 0.8 ml
Micro Bio-Spin P-6 Gel, SSC	7326200 7326201	25 x 0.8 ml 100 x 0.8 ml
Bio-Spin P-6 Gel, Tris	7326227	25 x 1.2 ml
Bio-Spin P-6 Gel, SSC	7326002 7326213	25 x 1.2 ml 1,000 x 1.2 ml
Micro Bio-Spin P-30 Gel, SSC	7326202 7326203 7326206	25 x 0.8 ml 100 x 0.8 ml 1,000 x 0.8 ml
Micro Bio-Spin P-30 Gel, Tris	7326223 7326226	25 x 0.8 ml 1,000 x 0.8 ml
Micro Bio-Spin P-30 Gel, Tris, RNase-free	7326250 7326251	25 x 0.8 ml 100 x 0.8 ml
Bio-Spin P-30 Gel, Tris	7326231 7326232	25 x 1.2 ml 100 x 1.2 ml
Bio-Spin P-30 Gel, SSC	7326006	100 x 1.2 ml
Mixed-Mode		
EconoFit CHT Type I, 40 µm	12009255 12009253 12009254	1 x 1 ml 1 x 5 ml 5 x 5 ml
Foresight CHT Type I, 40 µm	732-4735 732-4755	1 x 1 ml 1 x 5 ml
Foresight CHT Type I, 40 µm, RoboColumn Units	732-4822 732-4823	8 x 200 µl 8 x 600 µl
Foresight CHT Type I, 40 µm, Filter Plates	732-4716	2 x 96-well plates, 20 µl
EconoFit CHT Type II, 40 µm	12009259 12009257 12009258	1 x 1 ml 1 x 5 ml 5 x 5 ml
Foresight CHT Type II, 40 µm	732-4736 732-4756	1 x 1 ml 1 x 5 ml
Foresight CHT Type II, 40 µm, RoboColumn Units	732-4825 732-4826	8 x 200 µl 8 x 600 µl
Foresight CHT Type II, 40 µm, Filter Plates	732-4718	2 x 96-well plates, 20 µl
Bio-Scale CHT2-I	7510021	1 x 2 ml
Bio-Scale CHT5-I	7510023	1 x 5 ml
Bio-Scale CHT10-I	7510025	1 x 10 ml
Bio-Scale CHT20-I	7510027	1 x 20 ml

Description	Catalog #	Quantity
Prepacked Columns, continued		
EconoFit CHT Type I, 80 µm	12009256	1 x 1 ml
EconoFit CHT Type II, 80 µm	12009260	1 x 1 ml
EconoFit CHT XT, 40 µm	12009261	1 x 1 ml
Foresight CHT XT, 40 µm	12003150 12003149	1 x 1 ml 1 x 5 ml
EconoFit CFT Type II, 40 µm	12009252 12009240 12009251	1 x 1 ml 1 x 5 ml 5 x 5 ml
EconoFit MPC Type I, 40 µm	12009279	1 x 1 ml
Foresight MPC Type I, 40 µm	732-4737 732-4757	1 x 1 ml 1 x 5 ml
Foresight MPC Type I, 40 µm, RoboColumn Units	732-4828 732-4829	8 x 200 µl 8 x 600 µl
Foresight MPC Type I, 40 µm, Filter Plates	732-4785	2 x 96-well plates, 20 µl
EconoFit Nuvia aPrime 4A	12009280	1 x 1 ml
Foresight Nuvia aPrime 4A	12007392 12007393	1 x 1 ml 1 x 5 ml
Foresight Nuvia aPrime 4A RoboColumn Units	12007394 12007395	8 x 200 µl 8 x 600 µl
Foresight Nuvia aPrime 4A Filter Plates	12007411	2 x 96-well plates, 20 µl
EconoFit Nuvia cPrime	12009281	1 x 1 ml
Foresight Nuvia cPrime	732-4722 732-4742	1 x 1 ml 1 x 5 ml
Foresight Nuvia cPrime RoboColumn Units	732-4807 732-4808	8 x 200 µl 8 x 600 µl
Foresight Nuvia cPrime Filter Plates	732-4705	2 x 96-well plates, 20 µl
Ion Exchange		
ENrich Q	7800001 7800003	1 x 1 ml 1 x 8 ml
ENrich S	7800021 7800023	1 x 1 ml 1 x 8 ml
UNO Q1	7200001	1 x 1.3 ml
UNO Q6	7200003	1 x 6 ml
UNO Q12	7200005	1 x 12 ml
UNO Q Polishing	7200009	1 x 0.16 ml
UNO S1	7200021	1 x 1.3 ml
UNO S6	7200023	1 x 6 ml
UNO S12	7200025	1 x 12 ml
UNO S Polishing	7200029	1 x 0.16 ml
EconoFit Macro-Prep High Q	12009275 12009267 12009268 12009269	1 x 1 ml 5 x 1 ml 1 x 5 ml 5 x 5 ml
EconoFit Macro-Prep High Q-3HT	12009283	1 x 1 ml
EconoFit Macro-Prep DEAE	12009274 12009264 12009265 12009266	1 x 1 ml 5 x 1 ml 1 x 5 ml 5 x 5 ml
EconoFit Macro-Prep High S	12009276 12009270 12009271 12009272	1 x 1 ml 5 x 1 ml 1 x 5 ml 5 x 5 ml

Description	Catalog #	Quantity
Prepacked Columns, continued		
EconoFit Macro-Prep CM	12009273	1 x 1 ml
EconoFit UNOsphere Q	12009307	1 x 1 ml
	12009301	5 x 1 ml
	12009302	1 x 5 ml
	12009303	5 x 5 ml
Foresight UNOsphere Q	732-4732	1 x 1 ml
	732-4752	1 x 5 ml
Foresight UNOsphere Q RoboColumn Units	732-4819	8 x 200 µl
	732-4820	8 x 600 µl
Foresight UNOsphere Q Filter Plates	732-4714	2 x 96-well plates, 20 µl
EconoFit UNOsphere S	12009308	1 x 1 ml
	12009304	5 x 1 ml
	12009305	1 x 5 ml
	12009306	5 x 5 ml
Foresight UNOsphere S	732-4730	1 x 1 ml
	732-4750	1 x 5 ml
Foresight UNOsphere S RoboColumn Units	732-4813	8 x 200 µl
	732-4814	8 x 600 µl
Foresight UNOsphere S Filter Plates	732-4710	2 x 96-well plates, 20 µl
EconoFit Nuvia Q	12009290	1 x 1 ml
Foresight Nuvia Q	732-4721	1 x 1 ml
	732-4741	1 x 5 ml
Foresight Nuvia Q RoboColumn Units	732-4804	8 x 200 µl
	732-4805	8 x 600 µl
Foresight Nuvia Q Filter Plates	732-4703	2 x 96-well plates, 20 µl
EconoFit Nuvia HP-Q	12009282	1 x 1 ml
Foresight Nuvia HP-Q	12007020	1 x 1 ml
	12007021	1 x 5 ml
Foresight Nuvia HP-Q RoboColumn Units	12007013	8 x 200 µl
	12007014	8 x 600 µl
Foresight Nuvia HP-Q Filter Plates	12006908	2 x 96-well plates, 20 µl
EconoFit Nuvia S	12009291	1 x 1 ml
Foresight Nuvia S	732-4720	1 x 1 ml
	732-4740	1 x 5 ml
Foresight Nuvia S RoboColumn Units	732-4801	8 x 200 µl
	732-4802	8 x 600 µl
Foresight Nuvia S Filter Plates	732-4701	2 x 96-well plates, 20 µl
EconoFit Nuvia HR-S	12009284	1 x 1 ml
Foresight Nuvia HR-S	732-4723	1 x 1 ml
	732-4743	1 x 5 ml
Foresight Nuvia HR-S RoboColumn Units	732-4831	8 x 200 µl
	732-4832	8 x 600 µl
Foresight Nuvia HR-S Filter Plates	732-4707	2 x 96-well plates, 20 µl
Poly-Prep AG 1-X8, chloride, 100–200 mesh	7316211	50 x 2 ml

Description	Catalog #	Quantity
Prepacked Columns, continued		
Poly-Prep AG 1-X8, chloride, 200–400 mesh	7316212	50 x 2 ml
Poly-Prep AG 50W-X8, hydrogen, 100–200 mesh	7316213	50 x 2 ml
Poly-Prep AG 50W-X8, hydrogen, 200–400 mesh	7316214	50 x 2 ml
Poly-Prep AG 1-X8, formate, 200–400 mesh	7316221	50 x 2 ml
Hydrophobic Interaction		
EconoFit Macro-Prep t-Butyl	12009278	1 x 1 ml
EconoFit Macro-Prep Methyl	12009277	1 x 1 ml
Tagged Affinity		
EconoFit Nuvia IMAC	12009288	1 x 1 ml
	12009285	5 x 1 ml
	12009286	1 x 5 ml
	12009287	5 x 5 ml
Foresight Nuvia IMAC Ni-Charged	12004038	1 x 1 ml
	12004037	1 x 5 ml
Foresight Nuvia IMAC Ni-Charged RoboColumn Units	12004051	8 x 200 µl
	12004052	8 x 600 µl
Foresight Nuvia IMAC Ni-Charged Filter Plates	12004035	2 x 96-well plates, 20 µl
EconoFit Nuvia IMAC Uncharged	12009289	1 x 1 ml
EconoFit Profinity IMAC	12009298	5 x 1 ml
	12009299	1 x 5 ml
	12009300	5 x 5 ml
EconoFit Profinity GST	12009295	5 x 1 ml
	12009296	1 x 5 ml
	12009297	5 x 5 ml
EconoFit Profinity eXact	12009292	1 x 1 ml
	12009293	5 x 1 ml
	12009294	1 x 5 ml
Profinity eXact Mini Spin	1563007	10 x 0.1 ml
Immunoglobulin Affinity		
EconoFit UNOsphere SUPra	12009321	1 x 1 ml
	12009322	5 x 1 ml
	12009323	1 x 5 ml
EconoFit Affi-Prep Protein A	12009236	5 x 1 ml
	12009237	1 x 5 ml
Econo-Pac Protein A	7322022	5 x 2 ml
EconoFit Affi-Gel Blue	12009234	1 x 5 ml
	12009235	5 x 5 ml
EconoFit DEAE Affi-Gel Blue	12009262	1 x 5 ml
	12009263	5 x 5 ml

Ordering Information, continued

Description	Catalog #	Quantity
Empty Columns		
Econo Alpha Columns and Accessories		
Econo Alpha Column	12009463	6.6 x 50 mm
Econo Alpha Column	12009429	10 x 100 mm
Econo Alpha Column	12009430	15 x 100 mm
Econo Alpha Column	12009461	25 x 100 mm
Econo Alpha Column	12009462	10 x 500 mm
Econo Alpha Column	12009427	15 x 500 mm
Econo Alpha Column	12009428	25 x 500 mm
Econo Alpha Frits	12009464	15 mm, 5 µm pore size, pack of 20
Econo Alpha Frits	12009465	15 mm, 10 µm pore size, pack of 20
Econo Alpha Frits	12009466	15 mm, 20 µm pore size, pack of 20
Econo Alpha Frits	12009467	15 mm, 30 µm pore size, pack of 20
Econo Alpha Frits	12009468	15 mm, 50 µm pore size, pack of 20
Econo Alpha Frits	12009469	25 mm, 5 µm pore size, pack of 20
Econo Alpha Frits	12009470	25 mm, 10 µm pore size, pack of 20
Econo Alpha Frits	12009471	25 mm, 20 µm pore size, pack of 20
Econo Alpha Frits	12009472	25 mm, 30 µm pore size, pack of 20
Econo Alpha Frits	12009473	25 mm, 50 µm pore size, pack of 20
Econo Alpha Frits	12009474	10 mm, 5 µm pore size, pack of 20
Econo Alpha Frits	12009475	10 mm, 10 µm pore size, pack of 20
Econo Alpha Frits	12009476	10 mm, 20 µm pore size, pack of 20
Econo Alpha Frits	12009477	10 mm, 30 µm pore size, pack of 20
Econo Alpha Frits	12009478	10 mm, 50 µm pore size, pack of 20
Econo Alpha Frits	12009479	6.6 mm, 5 µm pore size, pack of 20
Econo Alpha Frits	12009480	6.6 mm, 10 µm pore size, pack of 20
Econo Alpha Frits	12009481	6.6 mm, 20 µm pore size, pack of 20
Econo Alpha Frits	12009482	6.6 mm, 30 µm pore size, pack of 20
Econo Alpha Frits	12009483	6.6 mm, 50 µm pore size, pack of 20
Econo Alpha Packing Sleeve	12009484	15 mm
Econo Alpha Packing Sleeve	12009485	25 mm
Econo Alpha Packing Sleeve	12009486	10 mm
Econo Alpha Packing Sleeve	12009487	6.6 mm
Econo Alpha O-Rings	12009488	15 mm

Description	Catalog #	Quantity
Empty Columns, continued		
Econo Alpha O-Rings	12009489	25 mm
Econo Alpha O-Rings	12009490	10 mm
Econo Alpha O-Rings	12009491	6.6 mm
Econo Alpha Adjustable Plunger	12009492	15 mm
Econo Alpha Adjustable Plunger	12009493	10 mm
Econo Alpha Adjustable Plunger	12009494	25 mm
Econo Alpha Column Extension	12009495	15 x 500 mm
Econo Alpha Column Extension	12009496	25 x 500 mm
Econo Alpha Column Extension	12009497	10 x 100 mm
Econo Alpha Column Extension	12009498	15 x 100 mm
Econo Alpha Column Extension	12009499	25 x 100 mm
Econo Alpha Column Extension	12009500	10 x 500 mm
Econo Alpha Column Extension	12009501	6.6 x 50 mm
Econo-Column Columns and Accessories		
Econo-Column Selection Pack A, 1 of each standard column	7376601	0.7 x 10, 20, 30 cm 1.5 x 30, 50 cm 2.5 x 20, 50 cm
Econo-Column Selection Pack B, 1 of each standard column	7376607	1.0 x 20, 30, 50 cm 1.5 x 20, 30, 50 cm
Econo-Column	7370507	0.5 x 5 cm, pack of 2
Econo-Column	7374506	0.5 x 5 cm, pack of 4
Econo-Column	7370512	0.5 x 10 cm, pack of 2
Econo-Column	7374516	0.5 x 15 cm, pack of 4
Econo-Column	7370517	0.5 x 15 cm, pack of 2
Econo-Column	7370522	0.5 x 20 cm, pack of 2
Econo-Column	7370707	0.7 x 5 cm, pack of 2
Econo-Column	7370712	0.7 x 10 cm, pack of 2
Econo-Column	7374711	0.7 x 10 cm, pack of 4
Econo-Column	7370717	0.7 x 15 cm, pack of 2
Econo-Column	7370732	0.7 x 30 cm, pack of 2
Econo-Column	7370752	0.7 x 50 cm, pack of 2
Econo-Column, Jacketed	7376108	0.7 x 15 cm, pack of 1
Econo-Column	7374006	1 x 5 cm, pack of 4
Econo-Column	7371007	1 x 5 cm, pack of 2
Econo-Column	7374011	1 x 10 cm, pack of 4
Econo-Column	7371012	1 x 10 cm, pack of 2
Econo-Column	7374021	1 x 20 cm, pack of 4
Econo-Column	7371022	1 x 20 cm, pack of 2
Econo-Column	7374031	1 x 30 cm, pack of 4
Econo-Column	7371032	1 x 30 cm, pack of 2
Econo-Column	7374051	1 x 50 cm, pack of 4
Econo-Column	7371052	1 x 50 cm, pack of 2
Econo-Column	7371091	1 x 100 cm, pack of 2
Econo-Column	7371093	1 x 120 cm, pack of 2
Econo-Column, Jacketed	7376116	1 x 15 cm, pack of 1
Econo-Column, Jacketed	7376131	1 x 30 cm, pack of 1
Econo-Column	7374150	1.5 x 5 cm, pack of 4
Econo-Column	7371507	1.5 x 5 cm, pack of 2

Description	Catalog #	Quantity
Empty Columns, continued		
Econo-Column	7374151	1.5 x 10 cm, pack of 4
Econo-Column	7371512	1.5 x 10 cm, pack of 2
Econo-Column	7374156	1.5 x 15 cm, pack of 4
Econo-Column	7371517	1.5 x 15 cm, pack of 2
Econo-Column	7374152	1.5 x 20 cm, pack of 4
Econo-Column	7371522	1.5 x 20 cm, pack of 2
Econo-Column	7374153	1.5 x 30 cm, pack of 4
Econo-Column	7371532	1.5 x 30 cm, pack of 2
Econo-Column	7374155	1.5 x 50 cm, pack of 4
Econo-Column	7371552	1.5 x 50 cm, pack of 2
Econo-Column	7371576	1.5 x 75 cm, pack of 2
Econo-Column	7371591	1.5 x 100 cm, pack of 2
Econo-Column	7371593	1.5 x 120 cm, pack of 2
Econo-Column	7371598	1.5 x 170 cm, pack of 2
Econo-Column, Jacketed	7376151	1.5 x 50 cm, pack of 1
Econo-Column	7372507	2.5 x 5 cm, pack of 2
Econo-Column	7374251	2.5 x 10 cm, pack of 4
Econo-Column	7372512	2.5 x 10 cm, pack of 2
Econo-Column	7374252	2.5 x 20 cm, pack of 4
Econo-Column	7372522	2.5 x 20 cm, pack of 2
Econo-Column	7374253	2.5 x 30 cm, pack of 4
Econo-Column	7372532	2.5 x 30 cm, pack of 2
Econo-Column	7372551	2.5 x 50 cm, pack of 2
Econo-Column	7372576	2.5 x 75 cm, pack of 2
Econo-Column	7372591	2.5 x 100 cm, pack of 2
Econo-Column	7372593	2.5 x 120 cm, pack of 2
Econo-Column	7375011	5 x 10 cm, pack of 1
Econo-Column	7375021	5 x 20 cm, pack of 1
Econo-Column	7375031	5 x 30 cm, pack of 1
Econo-Column	7375051	5 x 50 cm, pack of 1
Econo-Column	7375071	5 x 70 cm, pack of 1
Econo-Column Funnels	7310003	250 ml, pack of 5
Econo-Column Reservoir	7379112	500 ml
Econo-Column Reservoir	7379113	1 L
Flow Adaptor	7380014	1 cm inner diameter (ID), 1–7 cm length
Flow Adaptor	7380015	1 cm ID, 1–14 cm length
Flow Adaptor	7380016	1.5 cm ID, 1–14 cm length
Flow Adaptor	7380017	2.5 cm ID, 1–14 cm length
Flow Adaptor	7380018	5 cm ID, 1–14 cm length
Flow Adaptor Maintenance Kit	7380022	For 5 cm flow adaptor
	7380024	For 1 cm flow adaptor
	7380025	For 1.5 cm flow adaptor
	7380027	For 2.5 cm flow adaptor

Description	Catalog #	Quantity
Bio-Scale MT Columns and Accessories		
Bio-Scale MT2 Column	7510081	7 x 52 mm
Bio-Scale MT5 Column	7510083	10 x 64 mm
Bio-Scale MT10 Column	7510085	12 x 88 mm
Bio-Scale MT20 Column	7510087	15 x 113 mm
Replacement Parts Kit	7510091	For Bio-Scale MT2
	7510093	For Bio-Scale MT5
	7510095	For Bio-Scale MT10
	7510097	For Bio-Scale MT20
Bio-Spin, Mini Bio-Spin, and Micro Bio-Spin Columns		
Bio-Spin	7326008	100
Bio-Spin	7326025	1,000
Mini Bio-Spin	7326207	100
Micro Bio-Spin	7326204	100
End Caps for Bio-Spin, Micro Bio-Spin, Econo-Pac, and Poly-Prep	7311660	1,000
Econo-Pac Columns		
Econo-Pac	7321010	50
Econo-Pac	7321011	500
Poly-Prep Columns		
Poly-Prep	7311550	50
Poly-Prep	7311553	1,000
EconoFit Columns		
EconoFit	12009232	5 x 1 ml
EconoFit	12009233	5 x 5 ml

Visit [bio-rad.com/NGCsystem](https://www.bio-rad.com/NGCsystem) to learn more about the NGC Chromatography System.

Visit [bio-rad.com/MediaAndColumns](https://www.bio-rad.com/MediaAndColumns) for more information about Bio-Rad chromatography resins and columns.

BIO-RAD, AFFI-GEL, AFFI-PREP, AG, BIO-GEL, BIO-SPIN, CHELEX, CHT, ECONO-COLUMN, ECONO-PAC, MACRO-PREP, POLY-PREP, and UNO 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.

Profinity eXact vectors, tags, and resins are exclusively licensed under patent from Potomac Affinity Proteins. This product is intended for research purposes only. For any commercial use, commercial licenses are available by contacting Bio-Rad Laboratories, Inc., 6000 Alfred Nobel Drive, Hercules, CA 94547, attn.: Life Science Group Chromatography Marketing Manager, 1-800-4BIORAD (1-800-424-6723).

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



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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

