



## Resin and Column Selection Guide

The right column for the right resin

**BIO-RAD**

# Resin and Column Selection Guide

## Affinity | Desalting/Buffer Exchange



	Bio-Scale™ Mini Cartridges	Foresight™ Columns*	Bio-Scale MT Columns
<b>Main features</b>	Luer fitting with adaptors; easy self packing for automated or manual chromatography systems	Robust prepacked columns for high-throughput experiments	High pressure; acrylic-jacketed glass; adjustable
Maximum bed volume	1/5 ml	1/5 ml	2/5/10/20 ml
Bed height	0.56/1.3 cm	2/10 cm	5.2/6.4/8.8/11.3 cm
Maximum operating pressure	45 psi/0.3 MPa	435 psi/3 MPa	972 psi/6.7 MPa
Frit pore size	15–30 µm	7–12 µm	10 µm
Adaptors available for column	Yes	Yes	Yes
Available prepacked	Yes	Yes	Yes

### Chromatography Technique

Affinity Stationary Phase			
Nuvia™ IMAC	P	P	N
Profinity™ IMAC, Ni-Charged	P	–	N
Profinity IMAC	P	–	N
Profinity GST	P	–	N
Profinity eXact™	P	–	N
UNOsphere SUPrA™	P	P	N
Affi-Gel® Protein A	N	–	N
Affi-Prep® Protein A	P	–	R
DEAE Affi-Gel Blue	P	–	N
CM Affi-Gel Blue	R	–	N
Affi-Gel Blue	P	–	N
Affi-Gel Boronate	R	–	N
Affi-Prep Polymyxin	P	–	N
Affi-Gel Heparin	R	–	N
Affi-Gel HZ Hydrazide	R	–	N
Affi-Gel 102	R	–	N
Affi-Gel 10/15	R	–	N
Desalting/Buffer Exchange Stationary Phase			
Bio-Gel® P6 DG	P	–	N
Bio-Gel P30	R	–	N

\* Foresight Columns are available only in prepacked format.

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
<b>Main features</b>	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	N	N	N
Profinity IMAC, Ni-Charged	N	R	R
Profinity IMAC	N	R	R
Profinity GST	N	R	R
Profinity eXact	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	P	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 P6 DG	P	P	R
Bio-Gel P30	P	R	R

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

# Resin and Column Selection Guide

## Size Exclusion/Gel Filtration | Ion Exchange



	Bio-Scale™ Mini Cartridges	Foresight™ Columns*	Bio-Scale MT Columns
<b>Main features</b>	Luer fitting with adaptors; easy self packing for automated or manual chromatography systems	Robust prepacked columns for high-throughput experiments	High pressure; acrylic-jacketed glass; adjustable
Maximum bed volume	1/5 ml	1/5 ml	2/5/10/20 ml
Bed height	0.56/1.3 cm	2/10 cm	5.2/6.4/8.8/11.3 cm
Maximum operating pressure	45 psi/0.3 MPa	435 psi/3 MPa	971 psi/6.7 MPa
Frit pore size	15–30 µm	7–12 µm	10 µm
Adaptors available for column	Yes	Yes	Yes
Available prepacked	Yes	Yes	Yes

### Chromatography Technique

Ion Exchange Stationary Phase			
Macro-Prep® S	P	—	N
Macro-Prep Q	P	—	N
Macro-Prep Q-3HT	R	—	N
Macro-Prep 25 S	N	—	R
Macro-Prep 25 Q	N	—	R
Nuvia™ S	R	—	N
Nuvia Q	R	R	N
Nuvia HR-S	R	R	N
UNOsphere™ S	R	R	N
UNOsphere Q	P	R	N
AG®1	P	—	N
AG 2	N	—	N
AG 50W	N	—	N
AG 501	N	—	N
Bio-Rex™	N	—	N
Chelex®	N	—	N
ENrich™ S*,**	—	—	—
ENrich Q*,**	—	—	—
Bio-Rex 70	N	—	N
Bio-Rex RG 501-X8	N	—	N
Bio-Beads™ SM 2	N	R	N
Size Exclusion/Gel Filtration Stationary Phase			
Bio-Gel® P	P	—	N
Bio-Gel A	R	—	N
Bio-Beads S-X	R	—	O
ENrich 70*,**	—	—	—
ENrich 650*,**	—	—	—

\* Foresight and ENrich Columns are currently available only in prepacked format.

\*\* Ideal for use on an NGC Chromatography System or other medium pressure chromatography system.

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

# Resin and Column Selection Guide

## Size Exclusion/Gel Filtration | Ion Exchange



	Micro/Mini/Bio-Spin® Columns	Poly-Prep® and Econo-Pac® Columns	Glass Econo-Column® Columns
<b>Main features</b>	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
<b>Maximum bed volume</b>	0.8/1.2/1.6 ml	2/20 ml	1,374 ml
<b>Bed height</b>	2/2.8/3.7 cm	3.9/11.5 cm	5–170 cm
<b>Maximum operating pressure</b>	1,000 x g spin	14.7 psi/0.1 MPa	14.7 psi/0.1 MPa
<b>Frit pore size</b>	30 µm	30 µm	30 µm
<b>Adapters available for column</b>	N/A	Yes	Yes
<b>Available prepacked</b>	Yes	Yes	No

### Chromatography Technique

#### Ion Exchange Stationary Phase

Macro-Prep S	R	R	R
Macro-Prep Q	R	R	R
Macro-Prep Q-3HT	R	R	R
Macro-Prep 25 S	N	N	N
Macro-Prep 25 Q	N	N	N
Nuvia S	R	R	R
Nuvia Q	R	R	R
Nuvia HR-S	R	R	R
UNOsphere S	R	R	R
UNOsphere Q	R	R	R
AG 1	R	P	R
AG 2	R	R	R
AG 50W	R	P	R
AG 501	R	R	R
Bio-Rex	R	R	R
Chelex	R	R	R
ENrich S**	—	—	—
ENrich Q**	—	—	—
Bio-Rex 70	N	R	R
Bio-Rex RG 501-X8	N	R	R
Bio-Beads SM 2	N	N	N

#### 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
ENrich 70**	—	—	—
ENrich 650**	—	—	—

\* ENrich Columns are currently available only in prepacked format.

\*\* Ideal for use on an NGC Chromatography System or other medium pressure chromatography system.

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

# Resin and Column Selection Guide

## Hydrophobic Interaction | Mixed-Mode



	Bio-Scale™ Mini Cartridges	Foresight™ Columns*	Bio-Scale MT Columns
<b>Main features</b>	Luer fitting with adaptors; easy self packing for automated or manual chromatography systems	Robust prepacked columns for high-throughput experiments	High pressure; acrylic-jacketed glass; adjustable
<b>Maximum bed volume</b>	1/5 ml	1/5 ml	2/5/10/20 ml
<b>Bed height</b>	0.56/1.3 cm	N/A	5.2/6.4/8.8/11.3 cm
<b>Maximum operating pressure</b>	45 psi/0.3 MPa	435 psi/3 MPa	971 psi/6.7 MPa
<b>Frit pore size</b>	15–30 µm	7–12 µm	10 µm
<b>Adaptors available for column</b>	Yes	No	Yes
<b>Available prepacked</b>	Yes	Yes	Yes

### Chromatography Technique

#### Mixed-Mode Stationary Phase

CFT™, 40 µm	R	—	R
CFT II, 40 µm	R	—	R
CHT™ I, 20/40/80 µm	R	P	R
CHT II, 20/40/80 µm	R	P	R
MPC™	R	P	R
Bio-Gel® HT	N	N	N
Bio-Gel HTP	N	N	N
Nuvia™ cPrime™	R	P	N

#### Hydrophobic Interaction Stationary Phase

Macro-Prep® HIC (T-Butyl)	R	—	N
Bio-Beads® SM-2	R	—	N

\* Foresight Columns are available only in prepacked format.

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

# Resin and Column Selection Guide

## Hydrophobic Interaction | Mixed-Mode



	Micro/Mini/Bio-Spin® Columns	Poly-Prep® and Econo-Pac® Columns	Glass Econo-Column® Columns
<b>Main features</b>	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

Mixed-Mode Stationary Phase			
CFT, 40 µm	N	R	R
CFT II, 40 µm	N	R	R
CHT I, 20/40/80 µm	N	R	R
CHT II, 20/40/80 µm	N	R	R
MPC	N	R	R
Bio-Gel HT	N	R	R
Bio-Gel HTP	N	R	R
Nuvia cPrime	N	N	R
Hydrophobic Interaction Stationary Phase			
Macro-Prep HIC (T-Butyl)	R	R	R
Bio-Beads SM-2	R	R	R

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

## Ordering Information

### Prepacked Columns

Description	Catalog #	Quantity
<b>Gel Filtration/Desalting</b>		
ENrich SEC 70	7801070	24 ml
ENrich SEC 650	7801650	24 ml
	7324502	1 x 5 ml
	7324504	5 x 5 ml
Bio-Scale Mini	7325304	5 x 10 ml
Bio-Gel P-6	7325312	1 x 50 ml
	7325314	5 x 50 ml
<b>Mixed-Mode</b>		
Bio-Scale Mini CHT Type I	7324322	1 x 5 ml
	7324324	5 x 5 ml
Bio-Scale Mini CHT Type II	7324332	1 x 5 ml
	7324334	5 x 5 ml
Bio-Scale Mini CFT Type II	7324405	1 x 5 ml
	7324406	5 x 5 ml
	7324705	2 x 96-well plates
Foresight Nuvia cPrime	7324722	1 x 1 ml
	7324742	1 x 5 ml
<b>Ion Exchange</b>		
ENrich Q	7800001	1 ml
	7800003	8 ml
ENrich S	7800021	1 ml
	7800023	8 ml
UNO® Q	7200001	1 x 1.3 ml
	7200003	1 x 6 ml
	7200005	1 x 12 ml
UNO S	7200021	1 x 1.3 ml
	7200023	1 x 6 ml
	7200025	1 x 12 ml
Bio-Scale Mini Macro-Prep High Q	7324120	5 x 1 ml
	7324122	1 x 5 ml
	7324124	5 x 5 ml
Bio-Scale Mini Macro-Prep High S	7324130	5 x 1 ml
	7324132	1 x 5 ml
	7324134	5 x 5 ml
Bio-Scale Mini Macro-Prep DEAE	7324140	5 x 1 ml
	7324142	1 x 5 ml
	7324144	5 x 5 ml
Bio-Scale Mini UNOsphere S	7324110	5 x 1 ml
	7324112	1 x 5 ml
	7324114	5 x 5 ml
Foresight Nuvia Q	7324703	2 x 96-well plates
	7324721	1 x 1 ml
	7324741	1 x 5 ml
Foresight Nuvia S	7324720	1 x 1 ml
	7324740	1 x 5 ml
Bio-Scale Mini UNOsphere Rapid S	7324400	5 x 1 ml
	7324401	1 x 5 ml
	7324402	5 x 5 ml
Bio-Scale Mini UNOsphere Q	7324100	5 x 1 ml
	7324102	1 x 5 ml
	7324104	5 x 5 ml
<b>Tagged Affinity</b>		
Bio-Scale Mini Nuvia IMAC	7800811	1 x 5 ml
	7800812	5 x 5 ml
	7324620	5 x 1 ml
Bio-Scale Mini Profinity GST	7324622	1 x 5 ml
	7324624	5 x 5 ml
Bio-Scale Mini Profinity eXact	7324646	2 x 1 ml
	7324647	4 x 1 ml
	7324648	1 x 5 ml

### Prepacked Columns

Description	Catalog #	Quantity
<b>Immunoglobulin Affinity</b>		
Bio-Scale Mini UNOsphere SUPra Protein A	7324201	5 x 1 ml
	7324202	1 x 5 ml
Bio-Scale Mini Affi-Prep Protein A	7324600	5 x 1 ml
	7324602	1 x 5 ml
Bio-Scale Mini Affi-Gel Blue	7324642	1 x 5 ml
	7324644	5 x 5 ml

### Empty Columns

Description	Catalog #	Quantity
<b>Econo-Column Selection Packs</b>		
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
<b>Funnel and Reservoirs</b>		
Econo-Column Funnels	7310003	250 ml, 5/pk
Econo-Column Reservoir	7379112	500 ml
Econo-Column Reservoir	7379113	1 L
<b>Bio-Scale MT Columns</b>		
Bio-Scale MT2 Column	7510081	7 x 52 mm
Bio-Scale MT5 Column	7510083	10 x 64 mm
Bio-Scale MT20 Column	7510087	15 x 113 mm
Bio-Scale Mini Cartridge	7324660	1 ml, empty, 5/pk
Bio-Scale Mini Cartridge	7324661	5 ml, empty, 5/pk
<b>Bio-Spin, Mini Bio-Spin, and Micro Bio-Spin Columns</b>		
Bio-Spin Chromatography Columns	7326008	empty, 100
Bio-Spin Chromatography Columns	7326025	empty, 1,000
Micro Bio-Spin Chromatography Columns	7326204	empty, 100
Mini Bio-Spin Chromatography Columns	7326207	empty, 100
End Caps for Micro Bio-Spin Chromatography Columns	7311660	1,000
<b>Econo-Pac Columns</b>		
Econo-Pac Chromatography Columns	7321010	50
Econo-Pac Chromatography Columns	7321011	500
<b>Econo-Prep Columns</b>		
Poly-Prep Chromatography Columns	7311550	50
Poly-Prep Chromatography Columns	7311553	1,000

Visit [bio-rad.com/EmptyGlassColumns](http://bio-rad.com/EmptyGlassColumns) for a complete list of our empty glass Econo-Column Columns.

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

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

Profinity eXact vectors, tags, and resins, are licensed under patent rights from Potomac Affinity Proteins. For research use only. Any commercial use requires a license from Bio-Rad. 1-800-4BIORAD.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](http://bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 01 877 89019 Belgium 32 03 710 53 00 Brazil 55 11 3065 7550  
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 36 01 459 6192 Denmark 45 04 452 10 00 Finland 35 08 980 422 00  
France 33 01 479 593 00 Germany 49 089 3188 4393 Hong Kong 852 2789 3300 Hungary 36 01 459 6190 India 91 124 4029300  
Israel 972 03 963 6050 Italy 39 02 49486600 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 0 318 540 666  
New Zealand 64 9 415 2280 Norway 47 0 233 841 30 Poland 36 01 459 6191 Portugal 351 21 4727717 Russia 7 495 721 14 04  
Singapore 65 6415 3188 South Africa 36 01 459 6193 Spain 34 091 49 06 580 Sweden 46 08 555 127 00 Switzerland 41 0617 17 9555  
Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 01923 47 1301

