

---

# EconoFit Macro-Prep High Q, High S, DEAE, and CM Columns, 1 and 5 ml

## Instruction Manual

Catalog numbers

12009275  
12009267  
12009268  
12009269  
12009276  
12009270  
12009271  
12009272  
12009274  
12009264  
12009265  
12009266  
12009273

Please read the instructions in this manual prior to using EconoFit Macro-Prep High Q, High S, DEAE, and CM Columns. If you have any questions or require any further assistance, please contact your Bio-Rad Laboratories representative.

**BIO-RAD**

# Table of Contents

<b>Section 1</b>	Introduction . . . . .	1
<b>Section 2</b>	Product Information . . . . .	1
<b>Section 3</b>	Buffers and Methods . . . . .	2
<b>Section 4</b>	Preparing a Column and Subsequent Purification . . . . .	3
<b>Section 5</b>	Scaling Up . . . . .	4
<b>Section 6</b>	Regenerating, Cleaning, Sanitizing, and Storing Columns . . . . .	4
<b>Section 7</b>	Troubleshooting Guide . . . . .	5
<b>Section 8</b>	Ordering Information . . . . .	6
<b>Section 9</b>	Bibliography . . . . .	6

## Section 1

# Introduction

EconoFit Macro-Prep High Q, High S, DEAE, and CM Columns are convenient, disposable, prepacked low-pressure chromatography columns. EconoFit Columns offer both increased run-to-run uniformity and high purity of protein through the column design and novel resin technology. Compatible with aqueous buffers most commonly used for protein purification, EconoFit Columns offer improved performance for your protein separation needs.

EconoFit Macro-Prep High Q, High S, DEAE, and CM Columns are packed with Macro-Prep Ion Exchange Media. These media are based on hydrophilic, spherical, polymeric beads designed for the purification of proteins, nucleic acids, viruses, plasmids, and other macromolecules. Macro-Prep Beads are designed to provide medium capacity, low backpressure, and high productivity.

## Section 2

# Product Information

EconoFit Columns are disposable, easy-to-use, prepacked chromatographic columns supplied ready for use in convenient 1 and 5 ml sizes. They are quickly connected to liquid chromatography systems using 10-32 fittings. Columns are available for a variety of chromatographic techniques, including desalting (size exclusion [SEC]), ion exchange (IEX), affinity (AC), mixed-mode, and hydrophobic interaction chromatography (HIC). See Table 1 for specifications. Refer to [bio-rad.com/ResinsandColumns](http://bio-rad.com/ResinsandColumns) for a complete listing of items in the EconoFit Column product line.

**Table 1. EconoFit Macro-Prep High Q, High S, DEAE, and CM Column specifications.**

Property	Description
Size	1 and 5 ml bed volumes
Bed dimensions	1 ml: 25 mm length x 7 mm inner diameter 5 ml: 25 mm length x 16 mm inner diameter
Operational flow rates	1 ml: 1–6 ml/min (240–1,460 cm/hr) 5 ml: 5–15 ml/min (50–720 cm/hr)
Maximum operating pressure	72 psi
Fittings	10-32 (1/16"), female inlet and male outlet
Column material	Polypropylene
Frit material	High-density polyethylene
Shipping solution	20% ethanol
Storage conditions	20% ethanol
Autoclavability	Not autoclavable

Macro-Prep High Q, High S, DEAE, and CM Resins are also available in bottles. Refer to Ordering Information in section 8 of this manual. See Table 2 for specifications. Go to [bio-rad.com/ResinsandColumns](http://bio-rad.com/ResinsandColumns) for more information.

**Table 2. Macro-Prep High Q, High S, DEAE, and CM Resin specifications.**

Property	High Q	High S	DEAE	CM
Type of ion exchanger	Strong anion	Strong cation	Weak anion	Weak cation
Functional group	$-\text{N}^+(\text{CH}_3)_3$	$-\text{SO}_3^-$	$-\text{N}^+(\text{C}_2\text{H}_5)_2$	$-\text{COO}^-$
Nominal particle size	50 $\mu\text{m}$	50 $\mu\text{m}$	50 $\mu\text{m}$	50 $\mu\text{m}$
Nominal pore size	1,000 $\text{\AA}$	1,000 $\text{\AA}$	1,000 $\text{\AA}$	1,000 $\text{\AA}$
Total ionic capacity	400 $\pm$ 75 $\mu\text{eq/ml}$	160 $\pm$ 40 $\mu\text{eq/ml}$	175 $\pm$ 75 $\mu\text{eq/ml}$	210 $\pm$ 40 $\mu\text{eq/ml}$
Dynamic binding capacity*	$\geq$ 37 mg BSA/ml	$\geq$ 49 mg IgG*/ml	$\geq$ 30 mg BSA/ml	$\geq$ 25 mg hemoglobin/ml
Shipping counterion	$\text{Cl}^-$	$\text{Na}^+$	$\text{Cl}^-$	$\text{Na}^+$
Chemical stability				
1% SDS, 24 hr	Yes	Yes	Yes	Yes
6 M guanidine-HCl, 24 hr	Yes	Yes	Yes	Yes
pH stability	1–10	1–12	1–10	1–12
Antimicrobial agent	20% ethanol	20% ethanol	20% ethanol	20% ethanol
Regeneration	1–2 M NaCl	1–2 M NaCl	1–2 M NaCl	1–2 M NaCl
Storage conditions	20% ethanol	20% ethanol	20% ethanol	20% ethanol

\* 10% breakthrough capacity determined in a 1.1 x 20 cm column.

## Section 3

# Buffers and Methods

Ion exchange chromatography is usually performed using increasing salt gradients or pH gradients to elute the sample components. For best results, and increased column life, samples and buffers should be degassed and filtered through a 0.45  $\mu\text{m}$  filter.

Common buffers for cation and anion exchange chromatography are listed in Table 3.

An appropriate starting point for purifying samples is a linear gradient from 0–0.4 M NaCl spanning 1–20 column volumes at 120 cm/hr, 0.5 ml/min for the 1 ml column, and 2.5 ml/min for the 5 ml column. The separation can be optimized by changing the gradient profile. At the end of each run the column can be regenerated with 1.0 M NaCl followed by starting buffer. Return to the desired flow rate and proceed with the next separation.

**Table 3. Common buffers for ion exchange chromatography.**

Type of Buffering	
Cation	Ion Exchange Buffer Range, pH
Acetic acid	4.8–5.2
Citric acid	4.2–5.2
HEPES	7.6–8.2
Lactic acid	3.6–4.3
MES	5.5–6.7
MOPS	6.5–7.9
Phosphate	6.7–7.6
PIPES	6.1–7.5
Pivalic acid	4.7–5.4
TES	7.2–7.8
Tricine	7.8–8.9
Anion	Ion Exchange Buffer Range, pH
Bicine	7.6–9.0
Bis-Tris	5.8–7.2
Diethanolamine	8.4–8.8
Diethylamine	9.5–11.5
L-histidine	5.5–6.0
Imidazole	6.6–7.1
Pyridine	4.9–5.6
Tricine	7.8–8.9
Triethanolamine	7.3–8.0
Tris	7.5–8.0

## Section 4

# Preparing a Column and Subsequent Purification

EconoFit Macro-Prep High Q, High S, DEAE, and CM Columns contain 20% ethanol (v/v) as the storage solution. The fully hydrated support is ready to use after equilibrating the column in the buffer of choice. To perform buffer exchange, connect the column to a liquid chromatography system or peristaltic pump and condition it as instructed below:

1. Set pump flow rate to 3.0 ml/min (731 cm/hr) for the 1 ml column or 6.0 ml/min (288 cm/hr) for the 5 ml column.
2. Wash the column with degassed low salt buffer for 2 min.
3. Wash the column with degassed high salt buffer for 5 min.
4. Equilibrate the column with low salt buffer for 5 min.
5. Reduce the flow rate to the rate that will be used in the purification protocol.

## Sample Preparation

Proper pH and ionic strength are necessary for consistent and reproducible results. Sample can be exchanged into the starting buffer or diluted to the starting buffer concentration. This can be achieved by diluting the sample to the ionic strength of the starting buffer, dialyzing against the starting buffer, or exchanging it into the starting buffer. Buffer exchange can be accomplished using EconoFit Bio-Gel P6 Desalting Columns, Micro Bio-Spin P-6 or Micro Bio-Spin P-30 Columns, Bio-Spin P-6 or Bio-Spin P-30 Columns, Econo-Pac 10DG Desalting Columns, or Bio-Gel P-6DG Gel, as listed in Table 4. The choice of product will depend on the sample volume. All samples should be filtered through a 0.45  $\mu\text{m}$  filter prior to column application.

**Table 4. Product for buffer exchange.**

Sample Volume	Recommended Product	Use	Catalog #
10–75 $\mu\text{l}$	Micro Bio-Spin P-6 Column	Desalting proteins over 6 kD	7326221
10–75 $\mu\text{l}$	Micro Bio-Spin P-30 Column	Desalting proteins over 30 kD	7326223
50–100 $\mu\text{l}$	Bio-Spin P-6 Column	Desalting proteins over 6 kD	7326227
50–100 $\mu\text{l}$	Bio-Spin P-30 Column	Desalting proteins over 30 kD	7326231
100 $\mu\text{l}$ –3 ml	EconoFit Bio-Gel P6 Desalting Column	Desalting proteins over 6 kD	12009239
Up to 3 ml	Econo-Pac 10DG Desalting Columns	Desalting proteins over 6 kD	7322010
Unlimited	Bio-Gel P-6DG Gel	Desalting proteins over 6 kD	1500738

## Section 5 Scaling Up

EconoFit Columns are available in 1 and 5 ml formats. Macro-Prep High Q, High S, DEAE, and CM Resins are also available in various amounts, from 25 ml bottles to larger bulk quantities, for scaling up methods developed using the columns. For quick scale-up, two or three columns of the same type can be connected in series, so take care to maintain an overall system pressure  $\leq 72$  psi.

In addition, Bio-Rad carries an extensive line of empty chromatography columns from laboratory to process scale. Ask your local Bio-Rad representative or go to [bio-rad.com/ResinsandColumns](http://bio-rad.com/ResinsandColumns) for more information.

## Section 6 Regenerating, Cleaning, Sanitizing, and Storing Columns

### Regeneration

After each use, the column should be regenerated with the appropriate salt, in most cases 1–2 M NaCl in the presence of buffer. Wash with 2–4 column volumes of the buffered high salt solution. This reduces the potential for protein precipitation when selecting acid as a cleaning agent.

## Cleaning

After repeated use, an ion exchange column may require thorough cleaning and regeneration to remove bound contaminants. Acceptable cleaning-in-place (CIP) reagents include 1% acetic acid/1% phosphoric acid with 0.4 M NaCl, acetic acid (up to 30%), 1% Triton X-100, ethanol (up to 70%), isopropyl alcohol (up to 30%), 8 M urea, and 6 M guanidine-HCl. Any of these agents can be combined in an appropriate cleaning protocol. As a general guide, we recommend the following.

1. Use high salt buffer for regeneration, as above.
2. For aggregated or precipitated proteins, or when dirty feedstock (such as crude lysate) has been used, wash with 3–5 column volumes of 6 M guanidine-HCl or 8 M urea at 100 cm/hr.
3. For lipids or hydrophobically bound contaminants, wash with 0.1% Triton X-100 or 20–70% ethanol or isopropyl alcohol, or 1–30% acetic acid. Use 3–5 column volumes at 100 cm/hr.
4. Remove additional contaminants with 0.4 M NaCl in 1% acetic acid/1% phosphoric acid. Use 3–5 column volumes at 100 cm/hr.
5. If the column is to be used again immediately, wash with 2 column volumes of deionized water and 4–5 column volumes of starting buffer at 100 cm/hr. Check the conductivity and pH of the effluent to verify that the column is equilibrated in the starting buffer before loading the sample.

## Storage

After washing the columns with deionized water, EconoFit Ion Exchange Columns should be purged and stored with PBS containing 0.05% NaN<sub>3</sub>, or in 20% (v/v) ethanol solution, and capped for extended storage.

## Section 7 Troubleshooting Guide

Possible Causes	Possible Solutions
<b>Column Clogging or Slow Flow Rate</b>	
Particulates in sample	Filter all samples and buffers through 0.2 µm filter prior to application
<b>No Target Protein in Eluate</b>	
Low level of target	Check expression level of protein in starting SDS-PAGE material
<b>Precipitation during Purification</b>	
Binding capacity of column exceeded	Load less sample

## Section 8

# Ordering Information

Catalog #	Description
<b>EconoFit Macro-Prep High Q Columns</b>	
12009275	EconoFit Macro-Prep High Q Column, 1 x 1 ml column
12009267	EconoFit Macro-Prep High Q Columns, 5 x 1 ml columns
12009268	EconoFit Macro-Prep High Q Column, 1 x 5 ml column
12009269	EconoFit Macro-Prep High Q Columns, 5 x 5 ml columns
<b>EconoFit Macro-Prep High S Columns</b>	
12009276	EconoFit Macro-Prep High S Column, 1 x 1 ml column
12009270	EconoFit Macro-Prep High S Columns, 5 x 1 ml columns
12009271	EconoFit Macro-Prep High S Column, 1 x 5 ml column
12009272	EconoFit Macro-Prep High S Columns, 5 x 5 ml columns
<b>EconoFit Macro-Prep DEAE Columns</b>	
12009274	EconoFit Macro-Prep DEAE Column, 1 x 1 ml column
12009264	EconoFit Macro-Prep DEAE Columns, 5 x 1 ml columns
12009265	EconoFit Macro-Prep DEAE Column, 1 x 5 ml column
12009266	EconoFit Macro-Prep DEAE Columns, 5 x 5 ml columns
<b>EconoFit Macro-Prep CM Columns</b>	
12009273	EconoFit Macro-Prep CM Column, 1 x 1 ml column
<b>Macro-Prep High Q Resins</b>	
1580040	Macro-Prep High Q Support, 25 ml
1560040	Macro-Prep High Q Support, 100 ml
156-0041	Macro-Prep High Q Support, 500 ml
156-0042	Macro-Prep High Q Support, 5 L
156-0043	Macro-Prep High Q Support, 10 L
<b>Macro-Prep High S Resins</b>	
1580030	Macro-Prep High S Support, 25 ml
1560030	Macro-Prep High S Support, 100 ml
156-0031	Macro-Prep High S Support, 500 ml
156-0032	Macro-Prep High S Support, 5 L
156-0033	Macro-Prep High S Support, 10 L
<b>Macro-Prep DEAE Resins</b>	
1580020	Macro-Prep DEAE Support, 25 ml
1560020	Macro-Prep DEAE Support, 100 ml
156-0021	Macro-Prep DEAE Support, 500 ml
156-0022	Macro-Prep DEAE Support, 5 L
156-0023	Macro-Prep DEAE Support, 10 L
<b>Macro-Prep CM Resins</b>	
1580070	Macro-Prep CM Support, 25 ml
156-0070	Macro-Prep CM Support, 100 ml
156-0071	Macro-Prep CM Support, 500 ml
156-0073	Macro-Prep CM Support, 10 L

## Section 9

# Bibliography

- Gagnon P (1997). Avoiding instrument-associated aberrations in purification scale-up and scale-down. *BioPharm Int* 10, 42–45.
- Harris ELV and Angal S (1989). *Protein Purification Methods: A Practical Approach* (Oxford: IRL Press).
- Scopes RK (1987). *Protein Purification: Principles and Practice* (New York: Springer-Verlag).
- Snyder LR and Kirkland JJ (1979). *Introduction to Modern Liquid Chromatography* (New York: Wiley).



Bio-Rad, Bio-Gel, Bio-Spin, Econo-Pac, and Macro-Prep are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.  
All trademarks used herein are the property of their respective owner.



**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** 310 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 470 233 841 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

