



Bio-Scale DEAE Columns

Instruction Manual

**Catalog Number
751-0031, 751-0033,
751-0035, 751-0037**



Table of Contents

	Page
Section 1 Characteristics of the Bio-Scale DEAE2, DEAE5, DEAE10, and DEAE20 Anion Exchange Columns	1
1.1 Introduction	1
1.2 The DEAE Separation Medium	1
1.3 Connection to the BioLogic, FPLC® and HPLC Systems	1
Section 2 Use of the Bio-Scale DEAE Columns.....	2
2.1 Preparation for Initial Use	2
2.2 Buffer Selection.....	2
2.3 Sample Preparation	3
2.4 Elution Conditions.....	3
2.5 Use of Detergents	5
2.6 Chromatography Pre-Runs.....	5
2.7 Anomalous Chromatographic Behavior	5
Section 3 Care of the Bio-Scale DEAE Columns	6
3.1 Column Hygiene	6
3.2 Bed Height Adjustment.....	6
3.3 Frit Removal	6
3.4 Storage Conditions	6
3.5 References	8
Section 4 Product Information	8

Section 1

Characteristics of the Bio-Scale DEAE2, DEAE5, DEAE10, DEAE20 Anion Exchange Columns

1.1 Introduction

Bio-Scale prepacked ion exchange columns are designed to meet the needs of the bio-chromatographer for rapid and reproducible high resolution separations of biomolecules including proteins, peptides and polynucleotides. The availability of four column sizes (bed volumes of 2, 5, 10, and 20 ml) provides unrivaled flexibility for the economical and predictable scale up of separation and purification protocols without sacrificing resolution due to overloading.

1.2 The DEAE Separation Medium

Each column contains the spherical, rigid, and highly porous polymeric Macro-Prep[®] MP10 support derivatized with the weakly basic $-N^+(C_2H_5)_2$ diethylaminoethyl (DEAE) group. The 10 μ m particle size and narrow particle size distribution produces excellent resolution of biomolecules at high flow rates and with very low back-pressures.

The hydrophilic Macro-Prep DEAE support demonstrates extremely low non specific binding of biomolecules accompanied by high recovery of biological activity.

Stability of the MP10 DEAE Support

The columns are stable over the pH range 2-12, allowing for easy cleaning and regeneration. The MP10 support is compatible with aqueous solutions of guanidine-HCl and urea. Non-ionic and cationic detergents, and organic solvents such as methanol, ethanol and isopropanol, may also be used.

1.3 Connection to the BioLogic, FPLC[®] and HPLC Systems

Each Bio-Scale column is supplied with 0.02 inch ID (1/16 inch OD) Tefzel tubing, 1/4x28 flangeless fittings and end caps for connection to the BioLogic system. For connection to HPLC systems, use 10-32 fingertight fittings and M6 fittings for FPLC systems.

Instructions for assembling the fittings are included with the column.

Note: Bio-Rad Laboratories does not recommend nor warranty the use of the Bio-Scale columns with solvent delivery systems containing stainless steel parts when used with corrosive eluents containing eg halide salts. We recommend the use of inert, biocompatible (ceramic, PEEK, titanium) solvent delivery systems for maximum column life and recovery of sample biological activity.

Table 1. Column Characteristics

	DEAE2	DEAE5	DEAE10	DEAE20
Column volume (ml)	2	5	10	20
Recommended max. protein loading (mg/column)	20	50	100	200
Recommended flow rates (ml/min)	0.5 to 3.0	0.5 to 5.0	0.5 to 7.0	0.5 to 10.0
Static protein binding capacity (mg BSA/column)	50	125	250	500
Ionic capacity ($\mu\text{eg/ml}$)	200 ± 50	200 ± 50	200 ± 50	200 ± 50
Average particle size (μm)	10 ± 3.0	10 ± 3.0	10 ± 3.0	10 ± 3.0
Recommended buffer pH limits	5 - 9	5 - 9	5 - 9	5 - 9
Column dimensions (mm)	7 x 52	10 x 64	12 x 88	15 x 113
Maximum operating pressure (psi/bar)	1,000/69	750/50	600/40	500/34

Section 2 Use of the Bio-Scale DEAE Columns

2.1 Preparation for Initial Use

The columns are supplied in a storage buffer of 20% ethanol in water. The counter ion is OH⁻. Prior to initial use and after extended storage periods, each column should be conditioned as described below (steps 1-4). Always use HPLC grade reagents and filter and degas buffers). During this operation do not exceed more than **25%** of the maximum recommended flow-rates (see Table 1).

1. Wash with 5 column volumes of water.
2. Wash with 5 column volumes of 1M NaCl.
3. Wash with 5 column volumes of low ionic strength equilibration buffer. (*e.g.* 20 millimolar)
4. Wash with 5 column volumes of high ionic strength limit buffer. (*e.g.* 1 M NaCl)
5. Wash with 10 column volumes of low ionic strength equilibration buffer.

The column may now be further equilibrated in the starting buffer at the desired flow rate.

2.2 Buffer Selection

Table 2 lists commonly used buffers for anion exchange chromatography. These are cationic or zwitterionic buffers (it is important not to use anionic buffers which would interact with the anionic exchanger). The choice of whether to use an anion or cation exchanger is determined mainly by (a) the isoelectric point (pI) and, (b) the relationship between pH and the activity/stability of the protein of interest. Once the type of ion exchanger is determined, the choice of buffer and pH is also determined by the pH-activity relationship. As a general rule, the chosen buffer should be used in a pH range of $\text{pK}_a \pm 0.5$. This permits use of the lowest possible buffer concentration whilst maintaining maximum buffering capacity. **In any case, a minimum buffer concentration of 20 mM and pH range of 5 - 9 is recommended.**

As can be seen in Table 2, the pK_a and hence the pH of the buffer, changes with temperature. Therefore the pH of the buffer must be adjusted at the working temperature.

Table 2. Buffers for Anion-Exchange Chromatography

pH range	Buffer	Mwt	pK_a @25 °C	Counter-ion	ΔpK_a °C
5.0 - 6.0	Piperazine	86.1	5.7	Cl ⁻ /HCOO ⁻	-0.015
5.5 - 6.0	L-Histidine	155.2	6.15	Cl ⁻	
5.8 - 7.2	Bis-Tris	209.2	6.5	Cl ⁻	-0.017
6.4 - 7.3	Bis-Tris Propane	282.3	6.8, 9.0	Cl ⁻	
7.3 - 8.3	Triethanolamine	149.2	7.8	Cl ⁻ /CH ₃ COO ⁻	-0.020
7.6 - 8.6	Tris	121.1	8.1	Cl ⁻	-0.031
8.4 - 8.8	Diethanolamine	105.1	8.9	Cl ⁻	-0.025

Always use buffer components of the highest purity available as UV-absorbing impurities may cause baseline disturbances and interfere with the detection of protein peaks. Always filter and degas buffers.

2.3 Sample Preparation

Proper adjustment of the sample pH and ionic strength is critical for consistent and reproducible chromatography. For best results, the sample should be exchanged into the start buffer or diluted to the start buffer's concentration. Buffer exchange can be accomplished using Bio-Spin® 6 or Bio-Spin 30 columns, Econo-Pac® 10DG desalting columns, Bio-Gel® P-6DG size exclusion gel or the Econo-Pac P6 cartridge. The choice of product depends on sample volume. Always centrifuge or filter the sample (0.2-0.45 mm filter) to remove particulates. Application of turbid or lipid-containing samples may reduce the column lifetime. See Ref. 1 for further information about sample preparation.

Sample Load

The recommended sample load for each column is shown in Table 1. This amount may vary somewhat depending on the actual sample composition. We do not recommend overloading the column as both resolution and column lifetime will decrease. For larger loads, either change to a larger column or perform several chromatographic runs with a reduced loading. Ideally, samples should be bound in a concentrated zone at the top of the column. Higher sample loads produce a broad application zone in which components with less charge are displaced by more charged components. This may result in a shift of certain peaks to an earlier elution position in the gradient.

2.4 Elution Conditions

Separations by ion exchange are typically accomplished by increasing the salt concentration of the eluant either as a "step" or as a "continuous" gradient. For many separations, varying the pH of the elution buffer in addition to its salt concentration may be advantageous. Generally, it is best to choose initial pH and ionic strength conditions such that the protein of interest elutes early in the gradient. This is especially true for labile proteins or where a higher salt concentration is undesirable.

Choice of Elution Salt

Sodium or potassium chloride are the most common elution salts and are recommended for use with Bio-Scale DEAE columns. Other anions may be used and may show different selectivities based on their relative elution strengths.

These anions are shown below in increasing order of elution strengths:

Citrate>Sulfate>Iodide>Chloride>Formate>Acetate>Hydroxide.

See Ref. 2 for a more detailed explanation of ion selectivity.

Gradient Volumes and Salt Concentrations

As a starting point for developing a separation, we recommend using the Bio-Scale DEAE 2 column with a simple gradient profile over 48 ml.

Protocol: Use a flow rate of 2.0 ml/min. Following sample application, wash unbound proteins from the column with 6 ml (3 bed volumes) of Start buffer A. For elution, use a gradient volume of 20 ml (10 bed volumes) to a Cl^- concentration of 0.5 M (50% B). Follow this segment of the gradient by raising the salt concentration to 1.0 M (100% B) over 4 ml and then hold at 1.0 M for 6 ml before re-equilibrating the column with 12 ml of start buffer A. This gradient is shown schematically in Figure 1. Once an initial separation has been performed and the elution position of the protein of interest determined, the gradient composition and volume is adjusted to achieve maximum resolution. Normally, a gradient volume of 10 to 20 ml per ml of column bed volume is sufficient. The slope of the gradient will affect resolution. A steep gradient will result in relatively small peak volumes but short peak to peak distances. A shallower gradient normally gives greater resolution but peak volumes are larger.

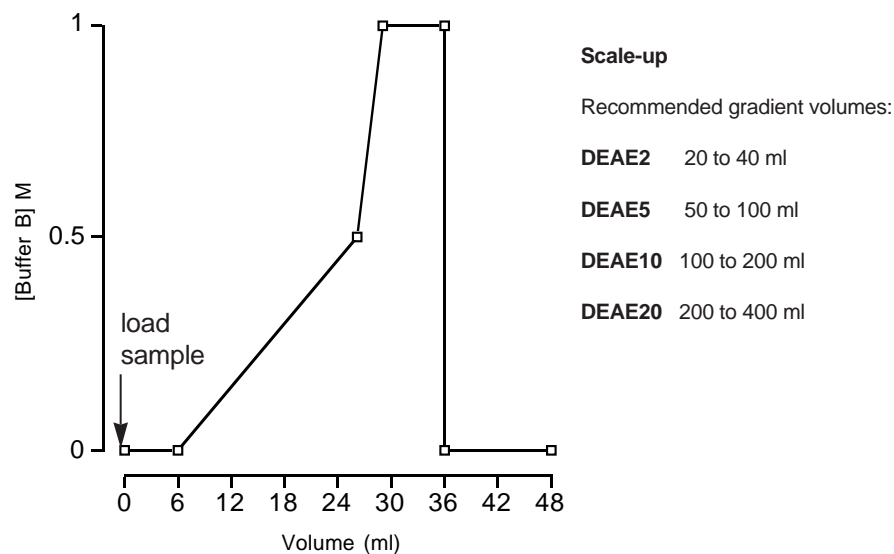


Fig. 1. Schematic gradient for separation on Bio-Scale DEAE2 column.

2.5 Use of Detergents

Cationic or non-ionic detergents may be used with the Bio-Scale DEAE support. We recommend the use of the reduced form of Triton[®] X-100 to minimize UV absorption artifacts. It is essential to thoroughly equilibrate the column with the detergent containing buffer prior to sample application. Pay particular attention to pH, which influences the solubility of the various classes of cationic detergents. Problems may arise when using salt gradient elution if the starting conditions include detergent below its critical micelle concentration (CMC). As the salt concentration (i.e. the counterion concentration) increases, the CMC drops and eventually micelles will form. This may cause a sudden increase in the UV baseline as the micelles themselves scatter light. We recommend using a concentration of detergent above the CMC during gradient elution. If subsequent chromatographic steps (eg. hydrophobic interaction) or other biochemical manipulations will be affected by the presence of the detergent, then the initial choice of detergent may be influenced by the availability of a protocol for its removal. See Refs. 1 and 3 for a more detailed explanation of the characteristics and use of detergents in biology and biochemistry and their removal from biological samples.

2.6 Chromatography Pre-Runs

Prior to loading the sample, we recommend running a blank gradient to check that the column is clean. In addition, detergents and other buffer components such as EDTA (at pH >2) and DTT (at pH > 8) will themselves elute as sharp, UV absorbing peaks during the gradient, complicating subsequent analysis of the chromatogram.

2.7 Anomalous Chromatographic Behavior

Certain proteins exhibit anomalous behavior on ion exchangers due to a non homogeneous charge distribution on their surface. Inclusion of zwitterions such as taurine and betaine (up to 200 mM) in the sample may improve the separation by delocalizing charges on such proteins. See Ref. 2 for further information.

Section 3 Care of the Bio-Scale DEAE Columns

3.1 Column Hygiene

Careful preparation (especially filtration) of the sample and the buffers will maintain the column performance and lifetime. Normally, washing with 1.0 M NaCl or KCl will remove most bound components. However, if there is a significant decrease in column performance (ie increasing backpressures or a significant drop in resolution) then a more extensive cleaning protocol as described below (steps 1-6) should be used. **Always reverse the flow during this procedure so tightly bound substances at the top of the column are quickly removed.**

During this operation do not exceed more than **25%** of the recommended maximum flow rates (see Table 1).

1. Wash with 2 Column Volumes (CV) of 1.0 M NaCl or KCl followed by 3 CV of water.
2. Wash with 2 x 1/2 CV of 0.5 M NaOH followed by 3 CV of water.
3. Wash with 2 x 1/2 CV of 2.0 M NaCl or KCl followed by 3 CV of water.
4. Wash with 1/2 CV of 30% Acetic Acid and 2 CV of 1M NaCl followed by 2 CV of water. Repeat.
5. If lipid contamination is a problem, wash with 1/2 CV of 50% methanol followed by 3 CV of water. Repeat.
6. Wash with 2 CV of 2.0 M NaCl or KCl or the salt containing the desired counter-ion.
7. Change the top frit, readjust the bed height to eliminate any voids and wash with 10 CV of equilibration buffer.

3.2 Bed Height Adjustment

Under certain conditions of buffer composition, high flow-rates or long term use, the resin bed may compress creating a void between the frit and the top of the bed. Normally, the void can be eliminated by turning the lock nut clockwise until the frit just touches the top of the bed. If the bed compresses at high flow rates, stop the pump and loosen the top fitting, then adjust the lock nut to remove the void, retighten the top fitting and resume pumping buffer.

Note: The Bio-Scale columns should not be run at their maximum flow rates for extended periods of time as this will cause excessive bed compression and high backpressures.

3.3 Frit Removal

The top frit and screen may need to be replaced after extensive column use or if increasing back pressures are noticed. Each column is supplied with a Frit Removal Tool, Polyethylene Frit and Distribution Screen. The screen assists in distributing the sample over the entire column surface and also acts as a pre filter for the frit. It should be replaced every time the frit is changed. Figure 2 describes in detail how to remove and change the top frit.

3.4 Storage Conditions

Prior to long term storage, the column should be cleaned as previously described and then washed with 3 column volumes of 20% ethanol. This will prevent microbial growth. Store the column in a safe place at room temperature. Never allow the column to freeze.

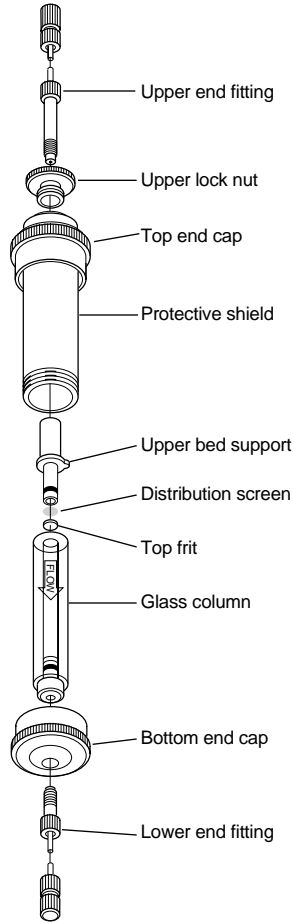


Fig. 2. Instructions for Removal of Top Frit.

1. Remove the upper and lower end fittings from the column.
2. Remove the upper lock nut from the top end cap.
3. Unscrew the bottom end cap from the protective shield and slide out the glass column.
4. Pull out the upper bed support from the glass column.
5. Using the frit removal tool, remove the distribution screen and frit by pressing the hook into the frit in a sideways motion with slight downward pressure.
6. Add 5 ml of buffer to the top of the resin bed. Place a new frit and screen into the tube. Insert the upper bed support and use it to carefully push down the frit and screen until it just touches the top of the resin bed.
7. Insert the glass column through the protective shield, into the top end-cap and align the bed support prong with the top end-cap guide. Attach the bottom end-cap to the shield. Insert the lock nut into the top end-cap and twist down until the plunger is at the top of the bed.
8. Attach column to a pump and pump buffer at 2.0 ml/min for 5 to 10 minutes. Inspect the column for voids and adjust accordingly.

3.5 References

1. Guide to Sample Preparation. Bio-Rad Laboratories Technical Bulletin # 1825.
2. W. Kopaciewicz, M. A. Rounds, J. Fausnaugh and F. E. Regnier (1983), Retention Model for High-performance Ion-Exchange Chromatography. *J. Chromatography* 266, 3-21.
3. A Guide to the Properties and Uses of Detergents in Biology and Biochemistry. J. Neugebauer (Calbiochem Biochemicals).

Section 4 Product Information

Catalog Number	Product Description
751-0031	Bio-Scale DEAE2 Column
751-0033	Bio-Scale DEAE5 Column
751-0035	Bio-Scale DEAE10 Column
751-0037	Bio-Scale DEAE20 Column
751-0039	Top-Off Resin Kit DEAE , includes 1 ml of MP10 DEAE resin, 5 frits and 5 distribution screens for each column diameter
751-0091	Bio-Scale 2 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover
751-0093	Bio-Scale 5 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover
751-0095	Bio-Scale 10 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover
751-0097	Bio-Scale 20 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover
751-0099	Bio-Scale Fittings Kit , includes 2 Super Flangeless Nuts (1/4 x 28 threads) and 4 ferrules, 2 Flangeless M6 nuts, 4 ferrules and 2 caps, 2 fingertight II fittings (10-32 threads)

Triton is a trademark of Rohm and Haas.

FPLC is a trademark of Pharmacia Biotech.

BIO-RAD**Bio-Rad
Laboratories****Life Science
Group**

2000 Alfred Nobel Drive
Hercules, California 94547
Telephone (510) 741-1000
Fax: (510) 741-5800

Australia, Bio-Rad Laboratories Pty Limited, Block Y Unit 1, Regents Park Industrial Estate, 391 Park Road, Regents Park, NSW 2143 • Phone 02-9414-2800 • Fax 02-9914-2888
Austria, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, 1130 Wien • Phone (1) 877 89 01 • Fax (1) 876 56 29
Belgium, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, 9810 Nazareth Eke • Phone 09-385 55 11 • Fax 09-385 65 54
Canada, Bio-Rad Laboratories (Canada) Ltd., 5671 McAdam Road, Mississauga, Ontario L4Z 1N9 • Phone (905) 712-2771 • Fax (905) 712-2990
China, Bio-Rad Laboratories, 14, Zhi Chun Road, Hai Dian District, Beijing 100088 • Phone (01) 2046622 • Fax (01) 2051876
Denmark, Bio-Rad Laboratories, Symbion Science Park, Fruebjergvej 3, DK-2100 Copenhagen • Phone 39 17 9947 • Fax 39 27 1698
Finland, Bio-Rad Laboratories, Business Center Länsikeskus, Pihatörmä 1A SF-02240, Espoo, • Phone 90 804 2200 • Fax 90 804 1100
France, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94 203 Ivry Sur Seine Cedex • Phone (1) 49 60 68 34 • Fax (1) 46 71 24 67
Germany, Bio-Rad Laboratories GmbH, Heidemannstraße 164, D-80939 München/Postfach 450133, D-80901 München • Phone 089 31884-0 • Fax 089 31884-100
India, Bio-Rad Laboratories, C-248 Defence Colony, New Delhi 110 024 • Phone 91-11-461-0103 • Fax 91-11-461-0765
Italy, Bio-Rad Laboratories S.r.l., Via Cellini, 18/A, 20090 Segrate Milano • Phone 02-21609 1 • Fax 02-21609-399
Japan, Nippon Bio-Rad Laboratories, 7-18, Higashi-Nippori 5-Chome, Arakawa-ku, Tokyo 116 • Phone 03-5811-6270 • Fax 03-5811-6272
The Netherlands, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 0318-540666 • Fax 0318-542216
New Zealand, Bio-Rad Laboratories Pty Ltd., P. O. Box 100-051, North Shore Mail Centre, Auckland 10 • Phone 09-443 3099 • Fax 09-443 3097
Pacific, Bio-Rad Laboratories, Unit 1111, 11/F., New Kowloon Plaza, 38, Tai Kok Tsui Road, Tai Kok Tsui, Kowloon, Hong Kong • Phone 7893300 • Fax 7891257
Singapore, Bio-Rad Laboratories (Singapore) Ltd., 221 Henderson Rd #05-19, Henderson Building, Singapore 0315 • Phone (65) 272-9877 • Fax (65) 273-4835
Spain, Bio-Rad Laboratories, S. A. Avda Valdelaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 85 • Fax (91) 661 96 98
Sweden, Bio-Rad Laboratories AB, Gärdsvägen 7D, Box 1276, S-171 24 Solna • Phone 46-(0)8-735 83 00 • Fax 46-(0)8-735 54 60
Switzerland, Bio-Rad Laboratories AG, Kanalstrasse 17, Postfach, CH-8152 Glattbrugg • Phone 01-809 55 55 • Fax 01-809 55 00
United Kingdom, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Free Phone 0800 181134 • Fax 01442 259118