
UNO™ Q&S
Continuous Bed
Ion Exchange Column
Instruction Manual

Catalog Numbers
720-0001, 720-0003, 720-0005,
720-0021, 720-0023 and 720-0025



UNO Anion and Cation Exchange Columns

Introduction

UNO 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 three column sizes (bed volumes of 1.3, 6, and 12 ml) provides unrivaled flexibility for the economical and predictable scale up of separation and purification protocols without sacrificing resolution due to overloading.

The UNO Separation Medium

Each column contains a “Continuous Bed” support derivatized with either strongly basic (-N⁺(CH₃)₃) or strongly acidic -SO₃⁻ sulfonic groups. The novel support produces excellent resolution of biomolecules at high flow rates and with low back-pressures. The hydrophilic support demonstrates extremely low non-specific binding of biomolecules accompanied by high recovery of biological activity.

Stability of the UNO Support

The columns are stable over the pH range 2–12, allowing for easy cleaning and regeneration. The support is compatible with aqueous solutions of guanidine-HCl and urea. Detergents and organic solvents such as 100% methanol, 100% ethanol, 100% isopropanol, and 30% ACN may also be used.

Connection to Chromatography Systems

Each UNO column is supplied with a fittings kit containing two pieces of 0.02 inch ID (1/16 inch OD) Tefzel tubing and two 1/4 x 28 flangeless fittings. Assemble one 1/4 x 28 fitting onto one end of the 0.02” tubing and insert each fitting into the top and bottom of the column. For connection to your chromatography system, attach the appropriate fitting onto the bare end of the Tefzel tubing.

Note: Bio-Rad Laboratories does not recommend nor warranty the use of the UNO columns with solvent delivery systems containing stainless steel parts when used with corrosive eluents containing *e.g.* 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. UNO Q Column Characteristics

	Q-1	Q-6	Q-12
Column bed volume (ml)	1.3	6	12
Recommended max. protein loading (mg)	20	80	160
Recommended flow-rates (ml/min)	0.5 to 5.0	0.5 to 8.0	0.5 to 10.0
Column Dimensions (mm)	7 x 35	12 x 53	15 x 68
Maximum operating pressure (psi/MPa/bar)	700/4.5/48	700/4.5/48	700/4.5/48

Table 2. UNO S Column Characteristics

	S-1	S-6	S-12
Column volume (ml)	1.3	6	12
Recommended max. protein loading (mg)	20	80	160
Recommended flow-rates (ml/min)	0.5 to 6.0	0.5 to 8.0	0.5 to 10.0
Column Dimensions (mm)	7 x 35	12 x 53	15 x 68
Maximum operating pressure (psi/MPa/bar)	700/4.5/48	700/4.5/48	700/4.5/48

Preparation for Initial Use

The columns are supplied in a storage buffer of 0.1 M NaCl + 20% ethanol. The counter ion for the Q column is Cl⁻ and Na⁺ for the S column. Prior to initial use and after extended storage periods, each column should be conditioned as described below (steps 1–4). Always use HPLC grade reagent, and filter and degas the buffers. During this operation do not exceed the following flow rates:

S-1 or Q-1	1 ml/min
S-6 or Q-6	2 ml/min
S-12 or Q-12	3 ml/min

1. Wash with 5 column volumes of water. Elevated backpressures may occur when washing with deionized water. Do not exceed 700 psi.
2. Wash with 5 column volumes of low ionic strength start buffer [*e.g.* 20 mM Tris-HCl (Q) or 20 mM Sodium Phosphate (S)].
3. Wash with 5 column volumes of high ionic strength elution buffer (*e.g.* Starting buffer + 1.0 M NaCl).
4. Wash with 5 column volumes of low ionic strength equilibration buffer [*e.g.* 20 mM Tris-HCl (Q) or 20 mM Sodium Phosphate (S)].

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

Buffer Selection

Table 2 and 3 lists commonly-used buffers for anion and cation-exchange chromatography. 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 pH used should be within ± 0.5 pH units of the pK_a of the chosen buffer. This permits use of the lowest possible buffer concentration while maintaining maximum buffering capacity. In any case, a buffer concentration of 20 mM is recommended.

As can be seen in Table 2 and 3, 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	Δ pK _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
9.0 - 9.9	Ethanolamine	61.1	9.5	Cl ⁻	-0.029
9.8 - 10.3	1,3-diamino-propane	74.1	10.5	Cl ⁻	-0.026

Table 3. Buffers for Cation-Exchange Chromatography

pH range	Buffer	Mwt	pK _a @25 °C	Counter-ion	Δ pK _a /°C
3.6 - 4.3	Lactic acid	90.1	3.8	Na ⁺	
4.2 - 5.2	Citric acid	192.1	3.1	Na ⁺	
5.5 - 6.7	MES	195.2	6.1	Na ⁺	-0.011
6.1 - 7.5	PIPES	302.4	6.8	Na ⁺	-0.009
6.5 - 7.9	MOPS	209.3	7.2	Na ⁺	-0.006
6.7 - 7.6	Phosphate	120 (Monobasic) 142 (Dibasic)	7.2	Na ⁺	-0.003
6.8 - 8.2	TES	229.2	7.4	Na ⁺	-0.020
6.8 - 8.2	HEPES	238.3	7.5	Na ⁺	-0.014
7.4 - 8.8	Tricine	179.2	8.1	Na ⁺	-0.021

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.

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.

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 switch to a larger column or perform several chromatographic runs with a reduced loading. Due to the fast flow properties of the UNO gel, multiple runs may be performed in a very short period of time. 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 highly charged components. This may result in a shift of certain peaks to an earlier elution position in the gradient.

Elution Conditions

Separations by ion-exchange chromatography are typically accomplished by increasing the salt concentration of the eluent 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 UNO columns. Other ions may be used and may show different selectivities based on their relative elution strengths and chaotropic nature.

The following ions are shown below in order of elution strengths:

Cations for UNO S Columns

Barium > Calcium > Magnesium > Potassium > Sodium > Lithium

Anions for UNO Q Columns

Citrate>Sulfate>Iodide>Chloride>Formate>Acetate

See Reference. 1 for a more detailed explanation of ion selectivity in chromatographic separations.

Gradient Volumes & Salt Concentrations

As a starting point for developing a separation, we recommend using the UNO S or UNO Q columns with a simple gradient profile over 10 column volumes.

Protocol: Use a flow-rate within the specified range in Table 1. Following sample application, wash unbound proteins from the column with 4 column volumes of Start buffer A. For elution, use a gradient volume of 10 column volumes to a NaCl 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 column volumes and then hold at 1.0 M for 4 column volumes before re-equilibrating the column with 5 column volumes 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.

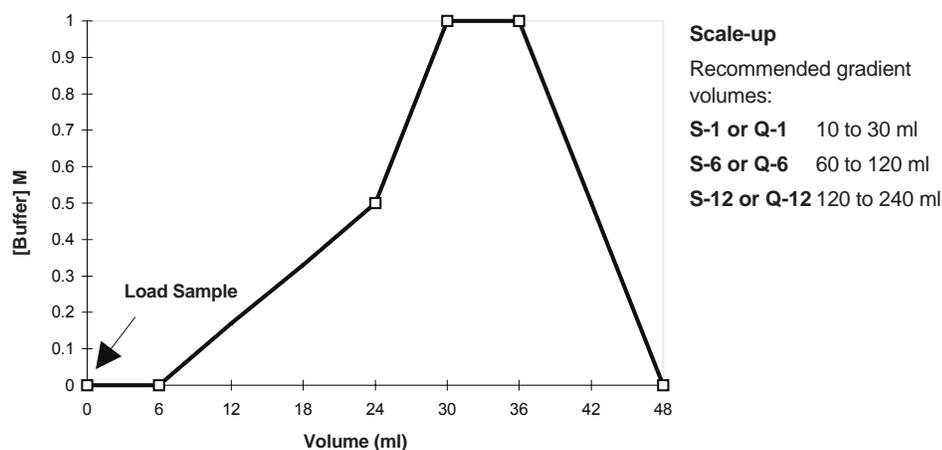


Fig. 1. Schematic gradient for separation on UNO S-1 or UNO Q-1 column.

Use of Detergents

Cationic or non-ionic detergents may be used with the UNO Q support. Anionic or non-ionic detergents may be used with the UNO S 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 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 (*e.g.* 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 Reference 2 for a more detailed explanation of the characteristics and use of detergents in biology and biochemistry and their removal from biological samples.

Chromatography Pre-Runs

Prior to loading the sample, it is recommended running a blank gradient to check that the column is clean. In addition, detergents and other buffer components will themselves elute as sharp, UV-absorbing peaks during the gradient, complicating subsequent analysis of the chromatogram.

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 (*i.e.* increasing backpressures or a significant drop in resolution) then a more extensive clean-

ing protocol as described below (steps 1–7) 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 the following maximum flow-rates.

S-1 or Q-1	1 ml/min
S-6 or Q-6	2 ml/min
S-12 or Q-12	3 ml/min

1. Wash with 2 column volumes of deionized water. Elevated backpressures may occur when washing with deionized water. Do not exceed 700 psi.
2. Wash with 2 column volumes of 2.0 M NaCl or KCl.
3. Wash with 1 column volume of 2.0 M NaOH followed by 3 column volumes of 0.1 M NaCl.
4. Wash with 1 column volume of 50% acetic acid followed by 3 column volumes of 0.1 M NaCl. Wash with 1 column volume of deionized water.
5. If lipid contamination is a problem, wash with 1 column volume of MeOH followed by 3 column volumes of 0.1 M NaCl.
6. Wash with 2 column volumes of 2.0 M NaCl or KCl or the salt containing the desired counter-ion.
7. Change the top frit and wash with 3 column volumes of start buffer.

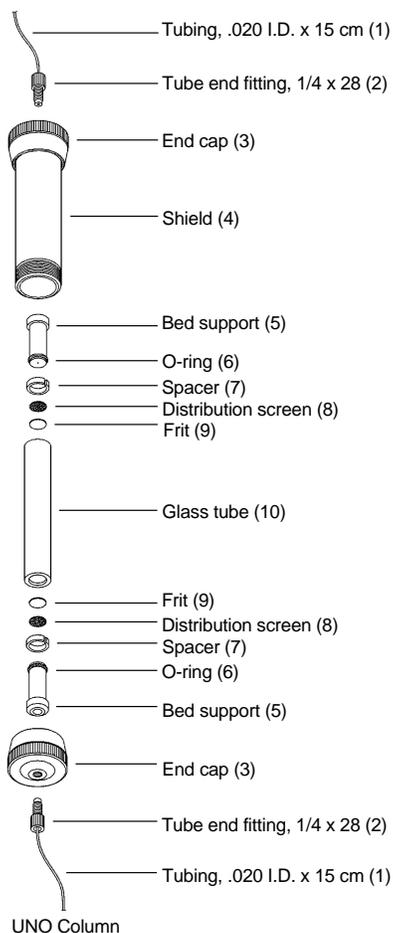


Fig. 1. Drawing of UNO column.

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 uniformly 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.

Column Repair

If the top of the resin bed becomes fouled and the above hygiene steps do not restore performance, then a few mm of the bed should be removed using the Bed Removal Tool. The tool will remove 4 mm of material. For detailed instructions on removing the bed material, see the instruction sheet included with the tool. Once fouled bed material is removed, insert a clean frit and distribution screen. **The dark brown ring on the top and bottom bed support must be removed before reinserting the bed supports into the glass tube. This procedure will compress the bed which is required to obtain proper performance.** This column repair procedure can only be performed once. If the back pressure of the column is still excessive, then the purchase of a UNO Replacement Column is recommended.

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.

References

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

Product Information

Catalog Number	Product Description
720-0001	UNO Q-1 Column
720-0003	UNO Q-6 Column
720-0005	UNO Q-12 Column
720-0009	UNO Q Polishing Column
720-0011	UNO Q-1 R Column , replacement column
720-0013	UNO Q-6 R Column , replacement column
720-0015	UNO Q-12 R Column , replacement column
720-0021	UNO S-1 Column
720-0023	UNO S-6 Column
720-0025	UNO S-12 Column

Catalog Number	Product Description
720-0029	UNO S Polishing Column
720-0031	UNO S-1 R Column , replacement column
720-0033	UNO S-6 R Column , replacement column
720-0035	UNO S-12 R Column , replacement column
751-0091	Bio-Scale 2 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover. Use this kit for the UNO S-1 or Q-1 Column.
751-0095	Bio-Scale 10 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover. Use this kit for the UNO S-6 or Q-6 Column.
751-0097	Bio-Scale 20 Replacement Part Kit , includes 5 frits, 5 distribution screens, 2 o-rings, 1 frit remover. Use this kit for the UNO S-12 or Q-12 Column.
751-0099	Bio-Scale Fittings Kit , includes 2 Super Flangeless Nuts (1/4 x 28 threads) and 6 ferrules, 2 Flangeless M6 Nuts, 4 ferrules and 2 caps, 2 Fingertight II fittings (10-32 threads). Use this kit for any UNO Column.
125-0561	Anion Exchange Standards , 6 pack
125-0562	Cation Exchange Standards , 6 pack

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, Pihatormä 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 31884-1000
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, Kanaistrasse 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