



## CHROMATOGRAPHY

# Nuvia™ Q Anion Exchange Resin

- Best-in-class binding capacity at fast flow rates
- Rapid mass transfer kinetics
- Efficient resolution of biomolecules from crude feedstreams
- Robust resin design providing competitive NaOH stability
- Full regulatory support

## High Binding Capacity and Resolution Capabilities for Downstream Processes

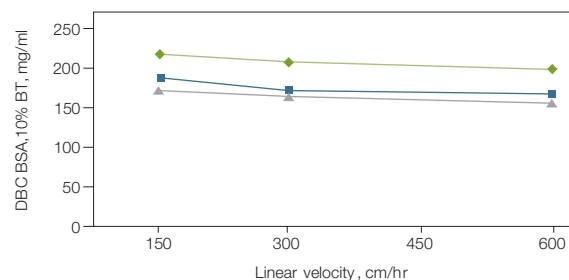
### Introduction

Nuvia Q is an ultra-high capacity, next-generation anion exchange resin built on the robust, industry-proven UNOsphere™ base matrix. Nuvia Q Resin delivers high binding capacity across a range of pH and flow rates, providing a wide experimental design space for process developers. This exceptional binding capacity makes it the anion exchanger of choice for biopharmaceuticals as upstream processes deliver higher concentrations of feed titers. Nuvia Q also delivers excellent performance for polishing applications. Nuvia Q can significantly improve productivity while contributing to reduced capital costs, space requirements, and cycle time for downstream biotherapeutic purification.

Nuvia Q is designed with proprietary surface extender technology on an optimized particle-size bead for improved flow and binding kinetics. The unique resin design gives Nuvia Q best-in-class dynamic binding capacity.

### High Capacity at Fast Flow Rate

Nuvia Q Resin is designed to meet the demanding pressure-flow requirements of downstream processes. With strong mechanical properties and optimized bead size distribution, Nuvia Q consistently delivers high binding capacity at fast linear flow velocities (Figure 1), allowing users the advantage of increased productivity without compromising binding capacity.



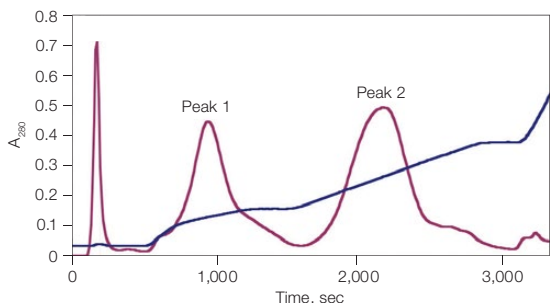
**Fig 1. Dynamic binding capacity vs. flow velocity of Nuvia Q Resin.** Each 1.1 cm column was packed to a 10.6 cm bed height with Nuvia Q (◆), agarose Q (■), or polymeric Q (▲) resin. Five mg/ml BSA in 20 mM Tris-HCl, pH 8.5, was loaded onto each column until 10% breakthrough was observed. BT, breakthrough; DBC BSA, dynamic binding capacity bovine serum albumin.

### Purification of Whey Proteins

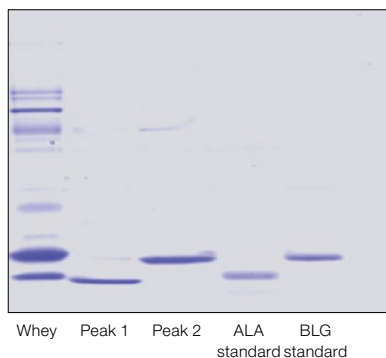
In addition to high dynamic binding capacity, Nuvia Q Resin also exhibits excellent baseline resolution of biomolecules with similar isoelectric points (pIs). Figure 2 shows the purification of  $\alpha$ -lactalbumin (ALA) and  $\beta$ -lactoglobulin (BLG), the major proteins of whey, accounting for approximately 75% of the total mass of crude whey.

The two major peaks were resolved using a segmented NaCl gradient when whey was used as a feedstream. SDS-PAGE analysis (Figure 3) of the fractions showed two protein peaks, one containing ALA (Peak 1) and the other containing BLG (Peak 2), with pIs of 4.7–5.1 and 5.2, respectively. Purities of the two purified proteins were estimated by densitometry to be at least 85%.





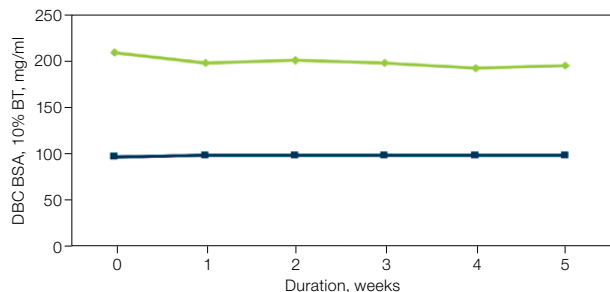
**Fig. 2. Purification of whey protein on Nuvia Q Resin.** Crude whey feed stock (8 mg) was loaded on a 7 x 27 mm column in 0.02 M sodium phosphate buffer, pH 6.0. The samples were eluted using a segmented gradient of 0–0.1 M NaCl, 0.1–0.3 M NaCl, and 0.3–0.9 M NaCl. Each fraction was 2 ml and analyzed by  $A_{280}$  (—); conductivity (—).



**Fig. 3. Analysis of whey purified on Nuvia Q Resin.** Fractions collected from the chromatography run shown in Figure 2 were separated by SDS-PAGE.

### Robust Performance

Nuvia Q Resin is produced by a validated manufacturing process to ensure strict batch-to-batch reproducibility. The chemical stability of Nuvia Q allows the resin to perform consistently with minimal change to dynamic binding capacity or recovery even after prolonged exposure to NaOH (Figure 4).



**Fig. 4. Stability of Nuvia Q Resin.** Results from base storage studies (in 1.0 N NaOH) show no loss in dynamic binding capacity and recovery. BSA binding capacity at 10% breakthrough (◆); recovery percentage (■). BT, breakthrough; DBC BSA, dynamic binding capacity bovine serum albumin.

### Properties of Nuvia Q Resin.

Property	Description
Type of ion exchanger	Strong anion
Functional group	$-N^+(CH_3)_3$
Total ionic capacity	100–170 $\mu\text{eq/ml}$
Dynamic binding capacity*	>170 mg/ml at 300 cm/hr
Shipping counterion	$\text{Cl}^-$
Particle size	$85 \pm 15 \mu\text{m}$
Recommended linear flow rate	50–600 cm/hr
Chemical stability	
1.0 N NaOH (20°C)	Up to 1 week
0.01 N NaOH (20°C)	Up to 5 years
Compression factor (settled bed volume/packed bed volume)	1.10–1.15
pH stability	
Short-term	2–14
Long-term	4–12
Shipping solution	20% ethanol + 1 M NaCl
Regeneration	1–2 M NaCl
Sanitization	0.5–1.0 N NaOH
Storage conditions	20% ethanol or 0.01 N NaOH

\* 10% breakthrough capacity determined with 5 mg/ml BSA in 20 mM Tris-HCl, pH 8.5.

### Technical Assistance

We have a dedicated applications team with extensive experience in process development that is available to answer questions and provide assistance in laboratory to scale-up design.

Regulatory support files are available upon request.

Bio-Rad Laboratories is an ISO 9001 registered corporation. For additional information and technical assistance, contact your local Bio-Rad office.

Visit [bio-rad.com/web/NuviaQ](http://bio-rad.com/web/NuviaQ) for more information about Bio-Rad's complete line of process chromatography resins.

### Ordering Information

Catalog #	Description
1560411	<b>Nuvia Q Resin</b> , 25 ml
1560413	<b>Nuvia Q Resin</b> , 100 ml
156-0415	<b>Nuvia Q Resin</b> , 500 ml
156-0417	<b>Nuvia Q Resin</b> , 10 L
732-4721	<b>Foresight™ Nuvia Q Column</b> , 1 ml
732-4741	<b>Foresight Nuvia Q Column</b> , 5 ml
732-4804	<b>Foresight Nuvia Q RoboColumn Unit</b> , 200 $\mu\text{l}$
732-4805	<b>Foresight Nuvia Q RoboColumn Unit</b> , 600 $\mu\text{l}$
732-4703	<b>Foresight Nuvia Q Plates</b> , 2 x 96-well, 20 $\mu\text{l}$

Larger volumes are available on request.

RoboColumn is a trademark of Atoll GmbH.



**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 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550  
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00  
France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 972 03 963 6050 Italy 39 02 216091 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 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04  
Singapore 65 6415 3188 South Africa 27 (0) 861 246 723 Spain 34 91 590 5200 Sweden 46 08 555 12700 Switzerland 41 026 674 55 05  
Taiwan 886 2 2578 7189 Thailand 66 662 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

