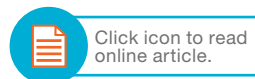


# Nuvia™ Chromatography Resins Publications List



Process Separations

Bulletin 6869



## Nuvia™ cPrime™ Hydrophobic Cation Exchange

### Wang R-Z et al. (2016).

New tetrapeptide ligands designed for antibody purification with biomimetic chromatography: Molecular simulation and experimental validation.  
Biochem Eng J 114, 191–201.



### Yan J et al. (2016).

Coadsorption of human immunoglobulin G and bovine serum albumin on a *p*-aminohippuric acid based mixed-mode resin.  
J Chem Eng Data 61, 151–159.



### Zhu M and Carta G (2016).

Protein adsorption equilibrium and kinetics in multimodal cation exchange resins.  
Adsorption 22, 165–179.



### Karkov HS et al. (2015).

Evaluation of selectivity in homologous multimodal chromatographic systems using in silico designed antibody fragment libraries.  
J Chromatogr A 1426, 102–109.



### Lončar N et al. (2015).

Mixed-mode resins: taking shortcut in downstream processing of raw-starch digesting  $\alpha$ -amylases.  
Sci Rep 5, 15772.



### Parimal S et al. (2015).

Interactions of multimodal ligands with proteins: insights into selectivity using molecular dynamics simulations.  
Langmuir 31, 7,512–7,523.



### Woo J et al. (2015).

Defining the property space for chromatographic ligands from a homologous series of mixed-mode ligands.  
J Chromatogr A 1407, 58–68.



### Woo J et al. (2015).

The effect of geometrical presentation of multimodal cation-exchange ligands on selective recognition of hydrophobic regions on protein surfaces.  
J Chromatogr A 1412, 33–42.



**BIO-RAD**



## Nuvia™ cPrime™ Hydrophobic Cation Exchange (continued)

### Srinivasan K et al. (2014).

Investigation into the molecular and thermodynamic basis of protein interactions in multimodal chromatography using functionalized nanoparticles.

Langmuir 30, 13,205–13,216.



### Yan J et al. (2014).

Protein adsorption behavior and immunoglobulin separation with a mixed-mode resin based on *p*-aminohippuric acid.

J Sep Sci 37, 2,474–2,480.



### Gao D et al. (2013).

Evaluating antibody monomer separation from associated aggregates using mixed-mode chromatography.

J Chromatogr A 1294, 70–75.



### Glaser V (2013).

Separation of therapeutic biomolecules: mixed-mode chromatography, layered bead designs, in silico modeling used at large scale.

Genetic Eng News 33, 33–34.



## Nuvia S Cation Exchange

### Guo J and Carta G (2015).

Unfolding and aggregation of monoclonal antibodies on cation exchange columns: effects of resin type, load buffer, and protein stability.

J Chromatogr A 1388, 184–194.



### Tao Y et al. (2014).

Evaluation of high-capacity cation exchange chromatography for direct capture of monoclonal antibodies from high-titer cell culture processes.

Biotechnol Bioeng 111, 1,354–1,364.



### Ng PK and Snyder MA (2012).

pH-based cation exchange chromatography in the capture and elution of monoclonal antibodies.

J Sep Sci 35, 29–35.



### Pérez Almodóvar EX et al. (2012).

Counterion effects on protein adsorption equilibrium and kinetics in polymer-grafted cation exchangers.

J Chromatogr A 1253, 83–93.



### Pérez Almodóvar EX et al. (2012).

Multicomponent adsorption of monoclonal antibodies on macroporous and polymer grafted cation exchangers.

J Chromatogr A 1264, 48–56.



### Pérez Almodóvar EX et al. (2011).

Protein adsorption and transport in cation exchangers with a rigid backbone matrix with and without polymeric surface extenders.

Biotechnol Prog 27, 1,264–1,272.





## Nuvia S Cation Exchange (continued)

**He X et al. (2010).**

Nuvia S Media.

Bioprocess Int 8, 59–61.



## Nuvia Q Anion Exchange

**Zhu M and Carta G (2014).**

Adsorption of polyethylene-glycolated bovine serum albumin on macroporous and polymer-grafted anion exchangers.

J Chromatogr A 1326, 29–38.



**Ng PK and Snyder MA (2013).**

Purification of  $\beta$ -lactoglobulin with a high-capacity anion exchanger: high-throughput process development and scale-up considerations.

J Sci Food Agric 93, 231–236.



**Nian R et al. (2013).**

Void exclusion of antibodies by grafted-ligand porous particle anion exchangers.

J Chromatogr A 1282, 127–132.



**Vetter TA et al. (2013).**

Mixed-beds of strong and weak anion exchange resins for protein separations with step-induced pH gradients.

Sep Sci Technol 49, 477–489.



**Ng PK and Snyder MA (2012).**

pH-based cation exchange chromatography in the capture and elution of monoclonal antibodies.

J Sep Sci 35, 29–35.



**Morrow KJ, Jr (2011).**

Strategies to advance mAb production: new approaches to process challenges help push the industry forward.

Genetic Eng News 31, 42–49.



**Pérez Almodóvar EX et al. (2011).**

Protein adsorption and transport in cation exchangers with a rigid backbone matrix with and without polymeric surface extenders.

Biotechnol Prog 27, 1,264–1,272.



Visit [bio-rad.com/web/ProcessPurification](http://bio-rad.com/web/ProcessPurification) to learn more about bioprocess production resins and their applications

**BIO-RAD**

**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 026674 55 05  
Taiwan 886 2 2578 7189 Thailand 66 662 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

