

# Macro-Prep<sup>®</sup> Chromatography Resins Publications List



Process Separations

Bulletin 6868



## Macro-Prep High Q Anion Exchange

### Mukhammadiev RS and Bagaeva TV (2015).

Isolation and structural characterization of *Rhizoctonia solani* fungal lectin.  
Res J Pharm Biol Chem Sci 6, 1,756–1,763.

### Bankston TE and Carta G (2010).

Apolipoprotein A-I(Milano) anion exchange chromatography: mass transfer and adsorption kinetics.  
Biotechnol J 5, 1,040–1,049.



### Liu B-X et al. (2008).

Purification and characterization of a leucine aminopeptidase from the skeletal muscle of common carp (*Cyprinus carpio*).  
Food Chem 108, 140–147.



### Abrunhosa L and Venâncio A (2007).

Isolation and purification of an enzyme hydrolyzing ochratoxin A from *Aspergillus niger*.  
Biotechnol Lett 29, 1,909–1,914.



### Thyer J et al. (2006).

Prion-removal capacity of chromatographic and ethanol precipitation steps used in the production of albumin and immunoglobulins.  
Vox Sang 91, 292–300.



### Song L (2006).

A soluble form of phosphatase in *Saccharomyces cerevisiae* capable of converting farnesyl diphosphate into E,E-farnesol.  
Appl Biochem Biotechnol 128, 149–158.



### Eon-Duval A and Burke G (2004).

Purification of pharmaceutical-grade plasmid DNA by anion-exchange chromatography in an RNase-free process.  
J Chromatogr B Analyt Technol Biomed Life Sci 804, 327–335.



### Song L (2003).

Detection of farnesyl diphosphate accumulation in yeast ERG9 mutants.  
Anal Biochem 317, 180–185.





## Macro-Prep High Q Anion Exchange (continued)

### Wilson MJ et al. (2001).

Removal of tightly bound endotoxin from biological products.  
J Biotechnol 88, 67–75.



## Macro-Prep DEAE Anion Exchange

### Liang TW et al. (2015).

Squid pen chitin chitoooligomers as food colorants absorbers.  
Mar Drugs 13, 681–696.



### Liang TW et al. (2015).

Chitinolytic bacteria-assisted conversion of squid pen and its effect on dyes and pigments adsorption.  
Mar Drugs 13, 4,576–4,593.



### Thinh PD et al. (2013).

Structural characteristics and anticancer activity of fucoidan from the brown alga *Sargassum mcclurei*.  
Mar Drugs 11, 1,456–1,476.



### Chanitnun K and Pinphanichakarn P (2012).

Glucose(xylose) isomerase production by *Streptomyces* sp. CH7 grown on agricultural residues.  
Braz J Microbiol 43, 1,084–1,093.



### Ermakova S et al. (2011).

Fucoidans from brown seaweeds *Sargassum hornery*, *Eclonia cava*, *Costaria costata*: structural characteristics and anticancer activity.  
Appl Biochem Biotechnol 164, 841–850.



### Sokolova RV et al. (2011).

Composition, structural characteristics, and antitumor properties of polysaccharides from the brown algae *Dictyopteris polypodioides* and *Sargassum* sp.  
Chemistry of Natural Compounds 47, 329.



### Wang et al. (2009).

Conversion and degradation of shellfish wastes by *Bacillus cereus* TKU018 fermentation for the production of chitosanases and bioactive materials.  
Biochem Eng J 48, 111–117.



### Chungool W et al. (2008).

Production, purification, and characterization of acetyl esterase from *Streptomyces* sp. PC22 and its action in cooperation with xylanolytic enzymes on xylan degradation.  
World J Microbiol Biotechnol 24, 549–556.





## Macro-Prep DEAE Anion Exchange (continued)

**Kim HJ et al. (2008).**

Membrane tetraheme cytochrome c(m552) of the ammonia-oxidizing *Nitrosomonas europaea*: a ubiquinone reductase.  
Biochemistry 47, 6,539–6,551.

**De Caro J et al. (2004).**

Characterization of pancreatic lipase-related protein 2 isolated from human pancreatic juice.  
Biochim Biophys Acta 1701, 89–99.

**Malmsten A et al. (2003).**

HIV-1 viral load determination based on reverse transcriptase activity recovered from human plasma.  
J Med Virol 71, 347–359.



## Macro-Prep High S Cation Exchange

**Leone S and Picone D (2016).**

Molecular dynamics driven design of pH-stabilized mutants of MNEI, a sweet protein.  
PLoS One 11, e0158372.

**Rodríguez-López A et al. (2016).**

Recombinant human N-acetylgalactosamine-6-sulfate sulfatase (GALNS) produced in the methylotrophic yeast *Pichia pastoris*.  
Sci Rep 6, 29329.

**Leone S et al. (2015).**

Acetate: friend or foe? Efficient production of a sweet protein in *Escherichia coli* BL21 using acetate as a carbon source.  
Microb Cell Fact 14, 106.

**Wang Y-J et al. (2015).**

Purification and enzymatic characterization of *E. coli* BL21 (DE3)/pET28a(+)-cr carbonyl reductase.  
Gao Xiao Hua Xue Gong Cheng Xue Bao/Journal of Chemical Engineering of Chinese Universities 29, 607–615.

**Davydov DR et al. (2013).**

CYP261 enzymes from deep sea bacteria: a clue to conformational heterogeneity in cytochromes P450.  
Biotechnol Appl Biochem 60, 30–40.

**Wang X et al. (2013).**

Preparation of the Mgm101 recombination protein by MBP-based tagging strategy.  
J Vis Exp 76, e50448.





## Macro-Prep High S Cation Exchange (continued)

**Golan-Mashiach M et al. (2012).**

Identification of CTCF as a master regulator of the clustered protocadherin genes.  
Nucleic Acids Res 40, 3,378–3,391.

**Lillehoj EP et al. (2012).**

NEU1 sialidase expressed in human airway epithelia regulates epidermal growth factor receptor (EGFR) and MUC1 protein signaling.  
J Biol Chem 287, 8,214–8,231.

**Park S and Lippard SJ (2012).**

Binding interaction of HMGB4 with cisplatin-modified DNA.  
Biochemistry 51, 6,728–6,737.

**Han L et al. (2008).**

A large form of secretogranin III functions as a sorting receptor for chromogranin A aggregates in PC12 cells.  
Mol Endocrinol 22, 1,935–1,949.

**Tsalkova TN et al. (2007).**

Mechanism of interactions of alpha-naphthoflavone with cytochrome P450 3A4 explored with an engineered enzyme bearing a fluorescent probe.  
Biochemistry 46, 106–119.



## Macro-Prep Hydrophobic Interaction Chromatography (HIC)

**Heger Z et al. (2015).**

Paramagnetic nanoparticles as a platform for FRET-based sarcosine picomolar detection.  
Sci Rep 5, 8868.

**Ribble W et al. (2015).**

Long-range PCR amplification of DNA by DNA polymerase III holoenzyme from *Thermus thermophilus*.  
Enzyme Res 2015, 837842.

**Wang Y-J et al. (2015).**

Purification and enzymatic characterization of *E. coli* BL21 (DE3)/pET28a(+)-cr carbonyl reductase.  
Gao Xiao Hua Xue Gong Cheng Xue Bao/Journal of Chemical Engineering of Chinese Universities 29, 607–615.

**Goto Y et al. (2011).**

KSAC, the first defined polyprotein vaccine candidate for visceral leishmaniasis.  
Clin Vaccine Immunol 18, 1,118–1,124.





## Macro-Prep Hydrophobic Interaction Chromatography (HIC) (continued)

### Michlmayr H et al. (2010).

Isolation and basic characterization of a beta-glucosidase from a strain of *Lactobacillus brevis* isolated from a malolactic starter culture. J Appl Microbiol 108, 550–559.



### Perumal SK et al. (2010).

Analysis of the DNA translocation and unwinding activities of T4 phage helicases. Methods 51, 277–288.



### Phogat S et al. (2008).

Analysis of the human immunodeficiency virus type 1 gp41 membrane proximal external region arrayed on hepatitis B surface antigen particles. Virology 373, 72–84.



### Soni B et al. (2008).

A novel method of single step hydrophobic interaction chromatography for the purification of phycocyanin from *Phormidium fragile* and its characterization for antioxidant property. Bioresour Technol 99, 188–194.



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

