

# Publication List: NGC Chromatography System Applications



Protein Purification

Bulletin 6843



Click icon to read  
online article.

## 2019

### Accumanno GM et al. (2019).

Purification and characterization of the thermostable protease produced by *Serratia grimesii* isolated from channel catfish.  
J Sci Food Agric 99, 2,428–2,437.



### Baruah GS et al. (2019).

Purification and characterization of phenoloxidase from the hemolymph of healthy and diseased *Antheraea assamensis* Helfer (*Lepidoptera: Saturniidae*): Effects of certain biological components and chemical agents on enzyme activity.  
Arch Insect Biochem Physiol 100, e21531.



### Becker W et al. (2019).

A fully automated three-step protein purification procedure for up to five samples using the NGC chromatography system.  
Protein Expr Purif 153, 1–6.



### de Araujo ED et al. (2019).

A functional in vitro assay for screening inhibitors of STAT5B phosphorylation.  
J Pharm Biomed Anal 162, 60–65.



### de la Cruz JJ et al. (2019).

Production of recombinant TSA-1 and evaluation of its potential for the immuno-therapeutic control of *Trypanosoma cruzi* infection in mice.  
Hum Vaccin Immunother 15, 210–219.



### Hoffmann D et al. (2019).

Downstream processing of Cry4AaCter-induced inclusion bodies containing insect-derived antimicrobial peptides produced in *Escherichia coli*.  
Protein Expr Purif 155, 120–129.



### Kern DM et al. (2019).

Cryo-EM structures of the DCPIB-inhibited volume-regulated anion channel LRRC8A in lipid nanodiscs.  
Elife 8, e42636.





## 2018

**Atykyan N et al. (2018).**

Isolation, purification, and investigation of some properties of glucose oxidase of the wood-degrading fungus *Lentinus (Panus) tigrinus* strain VKM F-3616D. BioResources 13 (issue 3).

**Baliban SM et al. (2018).**

Immunogenicity and induction of functional antibodies in rabbits immunized with a trivalent typhoid-invasive nontyphoidal *Salmonella* glycoconjugate formulation. Molecules 23, E1749.

**Bittel AM et al. (2018).**

Varied length Stokes shift BODIPY-based fluorophores for multicolor microscopy. Sci Rep 8, 4590.

**Chua LH et al. (2018).**

Process intensification of core streptavidin production through high-cell-density cultivation of recombinant *E. coli* and a temperature-based refolding method. J Biotechnol 276–277, 34–41.

**Dimond PF (2018).**

Scaleup issues remain front and center. Genetic Engineering & Biotechnology News 38, 16, 18.

**Frenzel E et al. (2018).**

In vivo selection of sfGFP variants with improved and reliable functionality in industrially important thermophilic bacteria. Biotechnol Biofuels 11, 8.

**Grzeszczuk MJ et al. (2018).**

Impact of selected amino acids of HP0377 (*Helicobacter pylori* thiol oxidoreductase) on its functioning as a CcmG (cytochrome c maturation) protein and Dsb (disulfide bond) isomerase. PLoS One 13, e0195358.

**Gumpena R et al. (2018).**

Crystal structure of the human dual specificity phosphatase 1 catalytic domain. Protein Sci 27, 561–567.

**Gumpena R et al. (2018).**

MBP-binding DARPins facilitate the crystallization of an MBP fusion protein. Acta Crystallogr F Struct Biol Commun 74, 549–557.

**Guo C et al. (2018).**

Molecular characterization of annexin B2, B3 and B12 in *Taenia multiceps*. Genes 9, E559.



**Krishnamurthy A and Belur PD (2018).**

A novel fibrinolytic serine metalloprotease from the marine *Serratia marcescens* subsp. *sakuensis*: Purification and characterization.  
Int J Biol Macromol 112, 110–118.

**Kudryakova IV et al. (2018).**

Structural and functional properties of antimicrobial protein L5 of *Lysobacter* sp. XL1.  
Appl Microbiol Biotechnol 102, 10,043–10,053.

**Kuo CY et al. (2018).**

Placental basement membrane proteins are required for effective cytotrophoblast invasion in a three-dimensional bioprinted placenta model.  
J Biomed Mater Res A 106, 1,476–1,487.

**Loi M et al. (2018).**

*In vitro* single and combined mycotoxins degradation by Ery4 laccase from *Pleurotus eryngii* and redox mediators.  
Food Control 90, 401–406.

**Luthra A et al. (2018).**

Structure and mechanism of a bacterial t6A biosynthesis system.  
Nucleic Acids Res 46, 1,395–1,411.

**Mathieu S et al. (2018).**

Ancient acquisition of “alginate utilization loci” by human gut microbiota.  
Sci Rep 8, 8075.

**Ou J et al. (2018).**

Bioprocess development of antibody-drug conjugate production for cancer treatment.  
PLoS One 13, e0206246.

**Pandey R et al. (2018).**

Differential diagnosis of otitis media with effusion using label-free Raman spectroscopy: A pilot study.  
J Biophotonics 11, e201700259.

**Rafikova O et al. (2018).**

Hemolysis-induced lung vascular leakage contributes to the development of pulmonary hypertension.  
Am J Respir Cell Mol Biol 59, 334–345.

**Retnoningrum DS et al. (2018).**

The first crystal structure of manganese superoxide dismutase from the genus *Staphylococcus*.  
Acta Crystallogr F Struct Biol Commun 74, 135–142.

**Selimoglu SM et al. (2018).**

Improved production of highly active and pure human creatine kinase MB.  
J Mol Microbiol Biotechnol 28, 28–36.



**Semenova MV et al. (2018).**

Cloning, isolation, and properties of a new homologous exoarabinase from the *Penicillium canescens* fungus.

Appl Biochem Microbiol 54, 387–395.

**Shi S et al. (2018).**

Human IGF-I propeptide A promotes articular chondrocyte biosynthesis and employs glycosylation-dependent heparin binding.

Biochim Biophys Acta Gen Subj 1862, 567–575.

**Smith RP et al. (2018).**

Production, biophysical characterization and initial crystallization studies of the N- and C-terminal domains of DsbD, an essential enzyme in *Neisseria meningitidis*.

Acta Crystallogr F Struct Biol Commun 74, 31–38.

**Toride KM and Brooks CL (2018).**

Epitope mapping of antibody-antigen interactions with X-ray crystallography.

Methods Mol Biol 1785, 13–27.

**Vafina G et al. (2018).**

Endonuclease from gram-negative bacteria *Serratia marcescens* is as effective as Pulmozyme in the hydrolysis of DNA in sputum.

Front Pharmacol 9, 114.

**Woznicka-Misaila A et al. (2018).**

Cell-free production, purification and characterization of human mitochondrial ADP/ATP carriers.

Protein Expr Purif 144, 46–54.

**Wu CH et al. (2018).**

Characterization of membrane-bound sulfane reductase: A missing link in the evolution of modern day respiratory complexes.

J Biol Chem 293, 16,687–16,696.

**Wu CH et al. (2018).**

Cytoplasmic and membrane-bound hydrogenases from *Pyrococcus furiosus*.

Methods Enzymol 613, 153–168.

**Wu CH et al. (2018).**

Improved production of the NiFe-hydrogenase from *Pyrococcus furiosus* by increased expression of maturation genes.

Protein Eng Des Sel 31, 337–344.

**Yang B et al. (2018).**

Comparative studies of the serum half-life extension of a protein via site-specific conjugation to a species-matched or -mismatched albumin.

Biomater Sci 6, 2,092–2,100.





## 2017

**Anderson GP et al. (2017).**

Thermal stabilization of anti- $\alpha$ -cobratoxin single domain antibodies.  
Toxicon 129, 68–73.

**Baena G et al. (2017).**

Salinity promotes opposite patterns of carbonylation and nitrosylation of  
C<sub>4</sub> phosphoenolpyruvate carboxylase in sorghum leaves.  
Planta 246, 1,203–1,214.

**Barth CW et al. (2017).**

Optimizing fresh specimen staining for rapid identification of tumor biomarkers  
during surgery.  
Theranostics 7, 4,722–4,734.

**Deguine J et al. (2017).**

Local TNFR1 signaling licenses murine neutrophils for increased TLR-dependent  
cytokine and eicosanoid production.  
J Immunol 198, 2,865–2,875.

**Fass R et al. (2017).**

Self-assembly of nanoscale particles with biosurfactants and membrane  
scaffold proteins.  
Biotechnol Rep 16, 1–4.

**Guo HF et al. (2017).**

A scalable lysyl hydroxylase 2 expression system and luciferase-based enzymatic  
activity assay.  
Arch Biochem Biophys 618, 45–51.

**Gutierrez PA et al. (2017).**

Differential effects of the dynein-regulatory factor Lissencephaly-1 on processive  
dynein-dynactin motility.  
J Biol Chem 292, 12,245–12,255.

**Jilek JL et al. (2017).**

Effects of microRNA-34a on the pharmacokinetics of cytochrome P450 probe drugs  
in mice.  
Drug Metab Dispos 45, 512–522.

**Liao JH et al. (2017).**

Acetylome of *Acinetobacter baumannii* SK17 reveals a highly-conserved modification  
of histone-like protein HU.  
Front Mol Biosci 4, 77.

**Maciejewska B et al. (2017).**

*Klebsiella* phages representing a novel clade of viruses with an unknown DNA  
modification and biotechnologically interesting enzymes.  
Appl Microbiol Biotechnol 101, 673–684.



**Maciejewska B et al. (2017).**

Modular endolysin of *Burkholderia* AP3 phage has the largest lysozyme-like catalytic subunit discovered to date and no catalytic aspartate residue.

Sci Rep 7, 14501.

**Mendes-Sousa AF et al. (2017).**

The sand fly salivary protein Lufaxin inhibits the early steps of the alternative pathway of complement by direct binding to the proconvertase C3b-B.

Front Immunol 8, 1065.

**Nilsson PH et al. (2017).**

Eculizumab-C5 complexes express a C5a neopeptide in vivo: Consequences for interpretation of patient complement analyses.

Mol Immunol 89, 111–114.

**Ogunmolu FE et al. (2017).**

Comparative insights into the saccharification potentials of a relatively unexplored but robust *Penicillium funiculosum* glycoside hydrolase 7 cellobiohydrolase.

Biotechnol Biofuels 10, 71.

**Rajiv C et al. (2017).**

The spliceosomal proteins PPIH and PRPF4 exhibit bi-partite binding.

Biochem J 474, 3,689–3,704.

**Retnoningrum DS et al. (2017).**

Structure-activity relationship of a recombinant hybrid manganese superoxide dismutase of *Staphylococcus saprophyticus*/*S. equorum*.

Int J Biol Macromol 98, 222–227.

**Selimoglu SM et al. (2017).**

Overcoming difficulties on synthesis of cardiac troponin-I.

Prep Biochem Biotechnol 47, 94–99.

**Sterky FH et al. (2017).**

Carbonic anhydrase-related protein CA10 is an evolutionarily conserved pan-neurexin ligand.

Proc Natl Acad Sci USA 114, E1253–E1262.

**Swendeman SL et al. (2017).**

An engineered S1P chaperone attenuates hypertension and ischemic injury.

Sci Signal 10, 2722.

**Urakami A et al. (2017).**

An envelope-modified tetra-valent dengue virus-like-particle vaccine has implications for Flavivirus vaccine design.

J Virol 91, e01181–e011817.

**Wang Y et al. (2017).**

A systematic protein refolding screen method using the DGR approach reveals that time and secondary TSA are essential variables.

Sci Rep 7, 9355.



**Wierzbicki IH et al. (2017).**

Functional and structural studies on the *Neisseria gonorrhoeae* GmhA, the first enzyme in the glycerol-mannose-heptose biosynthesis pathways, demonstrate a critical role in lipooligosaccharide synthesis and gonococcal viability.

Microbiologyopen 6, e00432.

**Wu R et al. (2017).**

Preparation of antioxidant peptides from salmon byproducts with bacterial extracellular proteases.

Mar Drugs 15, E4.

**Zhu Y et al. (2017).**

A cell-based, quantitative and isoform-specific assay for exchange proteins directly activated by cAMP.

Sci Rep 7, 6200.

**2016****Bonventre JA et al. (2016).**

Targeting an essential GTPase Opg for the development of broad-spectrum antibiotics.

PLoS One 11, e0148222.

**Cicéron F et al. (2016).**

Expression, purification and biochemical characterization of AtFUT1, a xyloglucan-specific fucosyltransferase from *Arabidopsis thaliana*.

Biochimie 128–129, 183–192.

**Çiftçi A et al. (2016).**

Development and validation of glycoprotein-based native-subunit vaccine for fish against *Aeromonas hydrophila*.

J Fish Dis 39, 981–992.

**Gong Z et al. (2016).**

Effect of a flexible linker on recombinant expression of cell-penetrating peptide fusion proteins and their translocation into fungal cells.

Mol Biotechnol 58, 838–849.

**Green AP et al. (2016).**

A chemically programmed proximal ligand enhances the catalytic properties of a heme enzyme.

J Am Chem Soc 138, 11,344–11,352.

**Holzer M et al. (2016).**

Refined purification strategy for reliable proteomic profiling of HDL<sub>2/3</sub>: Impact on proteomic complexity.

Sci Rep 6, 38533.



**Liebthal M et al. (2016).**

Redox-dependent conformational dynamics of decameric 2-cysteine peroxiredoxin and its interaction with cyclophilin 20-3.

Plant Cell Physiol 57, 1,415–1,425.

**Lunev S et al. (2016).**

Crystal structure of truncated aspartate transcarbamoylase from *Plasmodium falciparum*.

Acta Crystallogr F Struct Biol Commun 72, 523–533.

**Opoku-Temeng C and Sintim HO (2016).**

Inhibition of cyclic diadenylate cyclase, DisA, by polyphenols.

Sci Rep 6, 25445.

**Retnoningrum DS et al. (2016).**

Unique characteristics of recombinant hybrid manganese superoxide dismutase from *Staphylococcus equorum* and *S. saprophyticus*.

Protein J 35, 136–144.

**Tang W et al. (2016).**

High-level expression and characterization of the *Bacillus subtilis* subsp. *subtilis* str. BSP1 YwaD aminopeptidase in *Pichia pastoris*.

Protein Expr Purif 122, 23–30.

**Vafina G et al. (2016).**

A one-step protocol for chromatographic purification of non-recombinant exogenous bacterial enzyme: Nuclease of *Serratia marcescens*.

BioNanoScience 6, 335–337.

**Vdovin AS et al. (2016).**

Recombinant MHC tetramers for isolation of virus-specific CD8<sup>+</sup> cells from healthy donors: Potential approach for cell therapy of posttransplant cytomegalovirus infection.

Biochemistry 81, 1,371–1,383.

**Wang Z et al. (2016).**

Expression in *Pichia pastoris* and characterization of *Rhizomucor miehei* lipases containing a new propeptide region.

J Gen Appl Microbiol 62, 25–30.

**Williams PA et al. (2016).**

Phosphorylation of the RNA-binding protein Dazl by MAPKAP kinase 2 regulates spermatogenesis.

Mol Biol Cell 27, 2,341–2350.

**Zielke RA et al. (2016).**

Proteomics-driven antigen discovery for development of vaccines against gonorrhea.

Mol Cell Proteomics 15, 2,338–2,355.







## 2015

**An YD et al. (2015).**

Cloning, expression and purification of penicillin-binding protein 3 from *Pseudomonas aeruginosa* CMCC 10104. *Protein Expr Purif* 110, 37–42.

**Bocian-Ostrzycka KM et al. (2015).**

Functional and evolutionary analyses of *Helicobacter pylori* HP0231 (DsbK) protein with strong oxidative and chaperone activity characterized by a highly diverged dimerization domain. *Front Microbiol* 6, 1065.

**Cárdenas-Guerra RE et al. (2015).**

The recombinant prepro region of TvCP4 is an inhibitor of cathepsin L-like cysteine proteinases of *Trichomonas vaginalis* that inhibits trichomonal haemolysis. *Int J Biochem Cell Biol* 59, 73–83.

**Chen QX et al. (2015).**

A general approach to high-yield biosynthesis of chimeric RNAs bearing various types of functional small RNAs for broad applications. *Nucleic Acids Res* 43, 3,857–3,869.

**Degtyar E et al. (2015).**

Recombinant engineering of reversible cross-links into a resilient biopolymer. *Polymer* 69, 255–263.

**Frenkel ES and Ribbeck K (2015).**

Salivary mucins protect surfaces from colonization by cariogenic bacteria. *Appl Environ Microbiol* 81, 332–338.

**Ghodge SV and Raushel FM (2015).**

Discovery of a previously unrecognized ribonuclease from *Escherichia coli* that hydrolyzes 5'-phosphorylated fragments of RNA. *Biochemistry* 54, 2,911–2,918.

**Hernychová L et al. (2015).**

Recombinant expression, in vitro refolding and characterizing disulfide bonds of a mouse inhibitory C-type lectin-like receptor Nkrp1b. *Physiol Res* 64 (suppl 1), S85–S93.

**Kalia R et al. (2015).**

Structural and functional studies of membrane remodeling machines. *Methods Cell Biol* 128, 165–200.



**Lu J et al. (2015).**

One-step protein conjugation to upconversion nanoparticles.  
Anal Chem 87, 10,406–10,413.

**Ngo L et al. (2015).**

Effective quenchers are required to eliminate the interference of substrate:  
Cofactor binding in the HAT scintillation proximity assay.  
Assay Drug Dev Technol 13, 210–220.

**Padma Priya P et al. (2015).**

Characterization of precursor PfHsp60 in *Plasmodium falciparum* cytosol during its  
asexual development in human erythrocytes.  
PLoS One 10, e0136401.

**Roszczenko P et al. (2015).**

*Helicobacter pylori* HP0377, a member of the Dsb family, is an untypical multifunctional  
CcmG that cooperates with dimeric thioldisulfide oxidase HP0231.  
BMC Microbiol 15, 135.

**Su P et al. (2015).**

Isolation of Rhp-PSP, a member of YER057c/YjgF/UK114 protein family with antiviral  
properties, from the photosynthetic bacterium *Rhodospseudomonas palustris*  
strain JSC-3b.  
Sci Rep 5, 16121.

**Wang WP et al. (2015).**

Bioengineering novel chimeric microRNA-34a for prodrug cancer therapy: High-yield  
expression and purification, and structural and functional characterization.  
J Pharmacol Exp Ther 354, 131–141.

**Xiang DF et al. (2015).**

Interrogation of the substrate profile and catalytic properties of the phosphotriesterase  
from *Sphingobium* sp. strain TCM1: An enzyme capable of hydrolyzing  
organophosphate flame retardants and plasticizers.  
Biochemistry 54, 7,539–7,549.

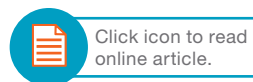
**2014****Huh I et al. (2014).**

In situ proteolysis, crystallization and preliminary X-ray diffraction analysis  
of a VHH that binds *Listeria* internalin B.  
Acta Crystallogr F Struct Biol Commun 70, 1,532–1,535.

**Li MM et al. (2014).**

Rapid production of novel pre-microRNA agent hsa-mir-27b in *Escherichia coli* using  
recombinant RNA technology for functional studies in mammalian cells.  
Drug Metab Dispos 42, 1,791–1,795.





---

**Sayeed SK et al. (2014).**

Identification of heat shock factor binding protein in *Plasmodium falciparum*.  
Malar J 13, 118.



---

**Traini M et al. (2014).**

Sphingomyelin phosphodiesterase acid-like 3A (SMPDL3A) is a novel nucleotide phosphodiesterase regulated by cholesterol in human macrophages.  
J Biol Chem 289, 32,895–32,913.



---

**Workman E et al. (2014).**

U1A regulates 3' processing of the survival motor neuron mRNA.  
J Biol Chem 289, 3,703–3,712.



---

Visit [bio-rad.com/web/ProteinPurification](http://bio-rad.com/web/ProteinPurification) for more information.



**Bio-Rad  
Laboratories, Inc.**

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.  
All trademarks used herein are the property of their respective owner.

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600  
**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 0 233 841 30  
**Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580  
**Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300  
**United Kingdom** 44 01923 47 1301

