



NGC Chromatography System Applications

Publications List

2021

Andreeva A et al. (2021).

Influence of carbohydrate additives on the growth rate of microalgae biomass with an increased carbohydrate content. *Mar Drugs* 19, 381.

Behar H et al. (2021).

Conservation of endo-glucanase 16 (EG16) activity across highly divergent plant lineages. *Biochem J* 478, 3,063–3,078.

Bergwik J et al. (2021).

Binding of the human antioxidation protein α 1-microglobulin (A1M) to heparin and heparan sulfate. Mapping of binding site, molecular and functional characterization, and co-localization *in vivo* and *in vitro*. *Redox Biol* 41, 101892.

Brouwer PJM et al. (2021).

Two-component spike nanoparticle vaccine protects macaques from SARS-CoV-2 infection. *Cell* 184, 1,188–1,200.e19.

Chan C et al. (2021).

The role of the *Moraxella catarrhalis* CopB protein in facilitating iron acquisition from human transferrin and lactoferrin. *Front Microbiol* 12, 714815.

Chatzileontiadou DSM et al. (2021).

Protein purification and crystallization of HLA-A*02:01 in complex with SARS-CoV-2 peptides. *Star Protoc* 2, 100635.

Deng L et al. (2021).

Bioengineered miR-124-3p prodrug selectively alters the proteome of human carcinoma cells to control multiple cellular components and lung metastasis *in vivo*. *Acta Pharm Sin B* 11, 3,950–3,965.

Felix J et al. (2021).

Structural and functional analysis of the *Francisella* lysine decarboxylase as a key actor in oxidative stress resistance. *Sci Rep* 11, 972.

Heymann D et al. (2021).

Structure of a consensus chitin-binding domain revealed by solution NMR.
J Struct Biol 213, 107725.

Kaku H et al. (2021).

Fas apoptosis inhibitory molecule blocks and dissolves pathological amyloid- β species.
Front Mol Neurosci 14, 750578.

Ke L et al. (2021).

Isolation and characterization of thermo-tolerant polyphenol oxidases in a black tea infusion.
Food Control 119, 107465.

Kern DM et al. (2021).

Cryo-EM structure of SARS-CoV-2 ORF3a in lipid nanodiscs.
Nat Struct Mol Biol 28, 573–582.

Krupka M et al. (2021).

Preparation and properties of recombinant *Clostridium ramosum* IgA proteinase. Isolation of Fc-SC and Fab fragments of human secretory IgA.
Protein Expr Purif 184, 105891.

Li PC et al. (2021).

In vivo fermentation production of humanized noncoding RNAs carrying payload miRNAs for targeted anticancer therapy.
Theranostics 11, 4,858–4,871.

Marcinkowski M et al. (2021).

Calmodulin as Ca²⁺-dependent interactor of FTO dioxygenase.
Int J Mol Sci 22, 10869.

Marivate A et al. (2021).

Recombinant expression, purification, and characterization of full-length human BST-2 from *Escherichia coli*.
Protein Expr Purif 188, 105969.

Martínez-Alarcón D et al. (2021).

Biochemical and structural studies of target lectin SapL1 from the emerging opportunistic microfungus *Scedosporium apiospermum*.
Sci Rep 11, 16109.

Messing S et al. (2021).

Production and membrane binding of N-terminally acetylated, C-terminally farnesylated and carboxymethylated KRAS4b.
Methods Mol Biol 2262, 105–116.

Ni K and Muegge K (2021).

LSH catalyzes ATP-driven exchange of histone variants macroH2A1 and macroH2A2.
Nucleic Acids Res 49, 8,024–8,036.

Parikh SJ et al. (2021).

Insights into the genetic variations of human cytochrome P450 2C9: Structural analysis, characterization and comparison. *Int J Mol Sci* 22, 10206.

Paxhia MD et al. (2021).

Functional characterization of the HMP-P synthase of *Legionella pneumophila* (Lpg1565). *Mol Microbiol* 115, 539–553.

Petrek H et al. (2021).

Single bioengineered ncRNA molecule for dual-targeting towards the control of non-small cell lung cancer patient-derived xenograft tumor growth.

Biochem Pharmacol 189, 114392.

Ren M et al. (2021).

Development of a super-sensitive diagnostic method for African swine fever using CRISPR techniques.

Virology 36, 220–230.

Ren M et al. (2021).

Early diagnosis of rabies virus infection by RPA-CRISPR techniques in a rat model.

Arch Virol 166, 1,083–1,092.

Si Y et al. (2021).

Anti-SSTR2 antibody-drug conjugate for neuroendocrine tumor therapy.

Cancer Gene Ther 28, 799–812.

2020

Ernst S et al. (2020).

Legionella effector AnkX displaces the switch II region for Rab1b phosphocolination.

Sci Adv 6, eaaz8041.

Fromell K et al. (2020).

Assessment of the role of C3(H₂O) in the alternative pathway.

Front Immunol 11, 530.

García-Cano I et al. (2020).

Purification and characterization of a phospholipid-hydrolyzing phosphoesterase produced by *Pediococcus acidilactici* isolated from Gouda cheese.

J Dairy Sci 103, 3,912–3,923.

Gulen B et al. (2020).

Identification of targets of AMPylating Fic enzymes by co-substrate-mediated covalent capture.

Nat Chem 12, 732–739.

Jeoung S et al. (2020).

Copper-binding energetics of amicyanin in different folding states.
Metallomics 12, 273–279.

Kaku H et al. (2020).

FAIM opposes aggregation of mutant SOD1 that typifies some forms of familial amyotrophic lateral sclerosis.
Front Neurosci 14, 110.

Ko MS et al. (2020).

Structurally plastic NEMO and oligomerization prone IKK2 subunits define the behavior of human IKK2:NEMO complexes in solution.

Biochim Biophys Acta Proteins Proteom 1868, 140526.

Li B et al. (2020).

Structural basis for pH gating of the two-pore domain K⁺ channel TASK2.

Nature 586, 457–462.

Nguyen TL et al. (2020).

The internal loops in the lower stem of primary microRNA transcripts facilitate single cleavage of human Microprocessor.
Nucleic Acids Res 48, 2,579–2,593.

Parikh SJ et al. (2020).

Structure of cytochrome P450 2C9*2 in complex with losartan: Insights into the effect of genetic polymorphism.

Mol Pharmacol 98, 529–539.

Placzkiewicz J et al. (2020).

Lactobacillus crispatus and its enolase and glutamine synthetase influence interactions between *Neisseria gonorrhoeae* and human epithelial cells.

J Microbiol 58, 405–414.

Sherekar M et al. (2020).

Biochemical and structural analyses reveal that the tumor suppressor neurofibromin (NF1) forms a high-affinity dimer.

J Biol Chem 295, 1,105–1,119.

Shibata K et al. (2020).

D-aspartate N-methyltransferase catalyzes biosynthesis of N-methyl-D-aspartate (NMDA), a well known selective agonist of the NMDA receptor, in mice.

Biochim Biophys Acta Proteins Proteom 1868, 140527.

Si Y et al. (2020).

Targeted exosomes for drug delivery: Biomanufacturing, surface tagging, and validation.

Biotechnol J 15, e1900163.

Tu MJ et al. (2020).

MicroRNA-1291-5p sensitizes pancreatic carcinoma cells to arginine deprivation and chemotherapy through the regulation of arginolysis and glycolysis.

Mol Pharmacol 98, 686–694.

Winters D et al. (2020).

Development of Bio-Rad NGC and GE ÄKTA pure systems for highly automated three column protein purification employing tandem affinity, buffer exchange and size exclusion chromatography.
Protein Expr Purif 165, 105497.

Yi W et al. (2020).

Bioengineered miR-328-3p modulates GLUT1-mediated glucose uptake and metabolism to exert synergistic antiproliferative effects with chemotherapeutics.
Acta Pharm Sin B 10, 159–170.

Yu Z et al. (2020).

Identification of protein-polysaccharide nanoparticles carrying hepatoprotective bioactives in freshwater clam (*Corbicula fluminea* Muller) soup.
Int J Biol Macromol 151, 781–786.

Zhu G et al. (2020).

TRIM11 prevents and reverses protein aggregation and rescues a mouse model of Parkinson's disease.
Cell Rep 33, 108418.

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.

Achek A et al. (2019).

Linear and rationally designed stapled peptides abrogate TLR4 pathway and relieve inflammatory symptoms in rheumatoid arthritis rat model.
J Med Chem 62, 6,495–6,511.

Ananthaswamy N et al. (2019).

A sequestered fusion peptide in the structure of an HIV-1 transmitted founder envelope trimer.
Nat Commun 10, 873.

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.

Fan X et al. (2019).

Non-affinity purification of a nanobody by void-exclusion anion exchange chromatography and multimodal weak cation exchange chromatography.
Sep Purif Technol 225, 88–96.

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.

Ilyina AP et al. (2019).

Analysis of serum albumin isoforms belonging to membranotropic homeostatic tissue-specific bioregulators isolated from different mammalian tissues.
Appl Biochem Microbiol 55, 355-359.

Kern DM et al. (2019).

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

Li X et al. (2019).

Bioengineered miR-27b-3p and miR-328-3p modulate drug metabolism and disposition via the regulation of target ADME gene expression.
Acta Pharm Sin B 9, 639–647.

Nikfarjam S et al. (2019).

Irreversible nature of mesoscopic aggregates in lysozyme solutions.
Colloid J 81, 546–554.

O'Neill LM et al. (2019).

Evaluation of protective and immune responses following vaccination with recombinant MIP and CPAF from *Chlamydia abortus* as novel vaccines for enzootic abortion of ewes.
Vaccine 37, 5,428–5,438.

Petrek H et al. (2019).

Bioengineering of a single long noncoding RNA molecule that carries multiple small RNAs.
Appl Microbiol Biotechnol 103, 6,107–6,117.

Wongpalee SP et al. (2019).

CryoEM structures of *Arabidopsis* DDR complexes involved in RNA-directed DNA methylation.
Nat Commun 10, 3,916.

Xu N et al. (2019).

Proteomics insight into the production of monoclonal antibody.
Biochem Eng J 145, 177–185.

Visit [bio-rad.com/NGC](https://www.bio-rad.com/NGC) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.
© 2022 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Website [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

