

# Publication List: NGC™ Chromatography System Applications



Protein Purification

Bulletin 6843



Click icon to read  
online article.

## 2016

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

Bonventre JA, Zielke RA, Korotkov KV, Sikora AE (2016). PLoS One 11, e0148222.



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

Çiftci A, Onuk EE, Çiftci G, Fındık A, Söğüt MÜ, Didinen BI, Aksoy A, Üstünakın K, Gülhan T, Balta F, Altun S (2016). J Fish Dis 39, 981–992.



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

Liebthal M, Strüve M, Li X, Hertle Y, Maynard D, Hellweg T, Viehhauser A, Dietz KJ (2016). Plant Cell Physiol [published online ahead of print February 12, 2016]. Accessed July 21, 2016.



### Inhibition of cyclic diadenylate cyclase, DisA, by polyphenols.

Opoku-Temeng C, Sintim HO (2016). Sci Rep 6, 25445.



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

Retnoningrum DS, Rahayu AP, Mulyanti D, Dita A, Valerius O, Ismaya WT (2016). Protein J 35, 136–144.



### Overcoming difficulties on synthesis of cardiac troponin-I.

Selimoglu SM, Kasap M, Akpınar G, Karadenizli A (2016). Prep Biochem Biotechnol [published online ahead of print April 12, 2016]. Accessed July 21, 2016.



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

Tang W, Li Z, Li C, Yu X, Wang F, Wan X, Wang Y, Ma L (2016). Protein Expr Purif 122, 23–30.





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

Wang Z, Lv P, Luo W, Yuan Z, He D (2016). *J Gen Appl Microbiol* 62, 25–30.



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

Zielke RA, Wierzbicki IH, Baarda BI, Gafken PR, Soge OO, Holmes KK, Jerse AE, Unemo M, Sikora AE (2016). *Mol Cell Proteomics* 15, 2,338–2,355.



**2015**

**Cloning, expression and purification of penicillin-binding protein 3 from *Pseudomonas aeruginosa* CMCC 10104.**

An YD, Du QZ, Tong LY, Yu ZW, Gong XW (2015). *Protein Expr Purif* 110, 37–42.



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

Bocian-Ostrzycka KM, Łasica AM, Dunin-Horkawicz S, Grzeszczuk MJ, Drabik K, Dobosz AM, Godlewska R, Nowak E, Collet JF, Jagusztyn-Krynicka EK (2015). *Front Microbiol* 6, 1065.



**The recombinant prepro region of TvCP4 is an inhibitor of cathepsin L-like cysteine proteinases of *Trichomonas vaginalis* that inhibits trichomonal haemolysis.**

Cárdenas-Guerra RE, Ortega-López J, Flores-Pucheta CI, Benítez-Cardoza CG, Arroyo R (2015). *Int J Biochem Cell Biol* 59, 73–83.



**A general approach to high-yield biosynthesis of chimeric RNAs bearing various types of functional small RNAs for broad applications.**

Chen QX, Wang WP, Zeng S, Urayama S, Yu AM (2015). *Nucleic Acids Res* 43, 3,857–3,869.



**Recombinant engineering of reversible cross-links into a resilient biopolymer.**

Degtyar E, Mlynarczyk B, Fratzl P, Harrington MJ (2015). *Polymer* 69, 255–263.

**Salivary mucins protect surfaces from colonization by cariogenic bacteria.**

Frenkel ES, Ribbeck K (2015). *Appl Environ Microbiol* 81, 332–338.



**Discovery of a previously unrecognized ribonuclease from *Escherichia coli* that hydrolyzes 5'-phosphorylated fragments of RNA.**

Ghodge SV, Raushel FM (2015). *Biochemistry* 54, 2,911–2,918.



**Recombinant expression, in vitro refolding and characterizing disulfide bonds of a mouse inhibitory C-type lectin-like receptor Nkrp1b.**

Hernychová L, Mrázek H, Ivanova L, Kukačka Z, Chmelík J, Novák P (2015). *Physiol Res* 64 Suppl 1, S85–93.



**Structural and functional studies of membrane remodeling machines.**

Kalia R, Talledge N, Frost A (2015). *Methods Cell Biol* 128, 165–200.

**One-step protein conjugation to upconversion nanoparticles.**

Lu J, Chen Y, Liu D, Ren W, Lu Y, Shi Y, Piper J, Paulsen I, Jin D (2015). *Anal Chem* 87, 10,406–10,413.

**Effective quenchers are required to eliminate the interference of substrate: cofactor binding in the HAT scintillation proximity assay.**

Ngo L, Wu J, Yang C, Zheng YG (2015). *Assay Drug Dev Technol* 13, 210–220.

**Characterization of precursor PfHsp60 in *Plasmodium falciparum* cytosol during its asexual development in human erythrocytes.**

Padma Priya P, Grover M, Tatu US, Natarajan V (2015). *PLoS One* 10, e0136401.

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

Roszczenko P, Grzeszczuk M, Kobierecka P, Wywiał E, Urbanowicz P, Wincek P, Nowak E, Jagusztyn-Krynicka EK (2015). *BMC Microbiol* 15, 135.

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

Su P, Feng T, Zhou X, Zhang S, Zhang Y, Cheng J, Luo Y, Peng J, Zhang Z, Lu X, Zhang D, Liu Y (2015). *Sci Rep* 5, 16121.

**Bioengineering novel chimeric microRNA-34a for prodrug cancer therapy: high-yield expression and purification, and structural and functional characterization.**

Wang WP, Ho PY, Chen QX, Addepalli B, Limbach PA, Li MM, Wu WJ, Jilek JL, Qiu JX, Zhang HJ, Li T, Wun T, White RD, Lam KS, Yu AM (2015). *J Pharmacol Exp Ther* 354, 131–141.

**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.**

Xiang DF, Bigley AN, Ren Z, Xue H, Hull KG, Romo D, Raushel FM (2015). *Biochemistry* 54, 7,539–7,549.

**2014****In situ proteolysis, crystallization and preliminary X-ray diffraction analysis of a VHH that binds listeria internalin B.**

Huh I, Gene R, Kumaran J, MacKenzie CR, Brooks CL (2014). *Acta Crystallogr F Struct Biol Commun* 70, 1,532–1,535.





**Rapid production of novel pre-microRNA agent hsa-mir-27b in *Escherichia coli* using recombinant RNA technology for functional studies in mammalian cells.**

Li MM, Wang WP, Wu WJ, Huang M, Yu AM (2014). Drug Metab Dispos 42, 1,791–1,795.



**Identification of heat shock factor binding protein in *Plasmodium falciparum*.**

Sayeed SK, Shah V, Chaubey S, Singh M, Alampalli SV, Tatu US (2014). Malar J 13, 118.



**Sphingomyelin phosphodiesterase acid-like 3A (SMPDL3A) is a novel nucleotide phosphodiesterase regulated by cholesterol in human macrophages.**

Traini M, Quinn CM, Sandoval C, Johansson E, Schroder K, Kockx M, Meikle PJ, Jessup W, Kritharides L (2014). J Biol Chem 289, 32,895–32,913.



**U1A regulates 3' processing of the survival motor neuron mRNA.**

Workman E, Veith A, Battle DJ (2014). 289, 3,703–3,712.



Visit [bio-rad.com/ProteinPurifications](http://bio-rad.com/ProteinPurifications) to learn more about the NGC Systems.



**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

