

Protein Sample Preparation Publications List



SureBeads™ Magnetic Beads System for Immunoprecipitation (IP)

Bulletin 7087



Click icon to read online article.

Selected Publications

Chen L and Brigstock D (2017).

Analysis of pathological activities of CCN proteins in fibrotic diseases: liver fibrosis. *Methods Mol Biol* 1489, 445–463.



Rathnasamy G et al. (2017).

Glutamate inhibits the pro-survival effects of insulin-like growth factor-1 on retinal ganglion cells in hypoxic neonatal rat retina. *Mol Neurobiol* 54, 3,453–3,464.



Anwar D et al. (2016).

p53 represses the transcription of *snRNA* genes by preventing the formation of little elongation complex. *BBA Gene Regul Mech* 1859, 975–982.



Azzariti A et al. (2016).

Hepatic stellate cells induce hepatocellular carcinoma cell resistance to sorafenib through the laminin-332/ α 3 integrin axis recovery of focal adhesion kinase ubiquitination. *Hepatology* 64, 2,103–2,117.



Böhm MR et al. (2016).

The pro-inflammatory role of high-mobility group box 1 protein (HMGB-1) in photoreceptors and retinal explants exposed to elevated pressure. *Lab Invest* 96, 409–427.



Czub B et al. (2016).

TOPORS, a dual E3 ubiquitin and SUMO1 ligase, interacts with 26S protease regulatory subunit 4, encoded by the PSMC1 gene. *PLoS One* 11, e0148678.



Elbahesh H et al. (2016).

Focal adhesion kinase (FAK) regulates polymerase activity of multiple influenza A virus subtypes. *Virology* 499, 369–374.



Fenstermaker RA et al. (2016).

Clinical study of a survivin long peptide vaccine (SurVaxM) in patients with recurrent malignant glioma. *Cancer Immunol Immunother* 65, 1,339–1,352.



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

An imbalance between specialized pro-resolving lipid mediators and pro-inflammatory leukotrienes promotes instability of atherosclerotic plaques.
Nat Commun 7, 12859.

**Hashimoto K et al. (2016).**

TAK1 regulates the Nrf2 antioxidant system through modulating p62/SQSTM1.
Antioxid Redox Signal 25, 953–964.

**Hoppstädter J et al. (2016).**

Induction of glucocorticoid-induced leucine zipper (GILZ) contributes to anti-inflammatory effects of the natural product curcumin in macrophages.
J Biol Chem 291, 22,949–22,960.

**Kim GT et al. (2016).**

Torilis japonica extract-generated intracellular ROS induces apoptosis by reducing the mitochondrial membrane potential via regulation of the AMPK-p38 MAPK signaling pathway in HCT116 colon cancer.
Intl J Oncol 49, 1,088–1,098.

**Kumari R et al. (2016).**

Movement protein of *cucumber mosaic virus* associates with apoplasmic ascorbate oxidase.
PLoS One 11, e0163320.

**Masatani T et al. (2016).**

Contribution of the interaction between the rabies virus P protein and I-kappa B kinase ϵ to the inhibition of type I IFN induction signalling.
J Gen Virol 97, 316–326.

**Menden HL et al. (2016).**

NADPH oxidase 2 regulates LPS-induced inflammation and alveolar remodeling in the developing lung.
Am J Respir Cell Mol Biol 55, 767–778.

**Muraki M (2016).**

Preparation of a functional fluorescent human Fas ligand extracellular domain derivative using a three-dimensional structure guided site-specific fluorochrome conjugation.
SpringerPlus 5, 997.

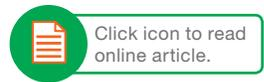
**Nowotarski SL et al. (2016).**

Destabilization of the ornithine decarboxylase mRNA transcript by the RNA-binding protein tristetraprolin.
Amino Acids 48, 2,302–2,311.

**Ozaki H et al. (2016).**

Ca²⁺/calmodulin-dependent protein kinase phosphatase (CaMKP/PPM1F) interacts with neurofilament L and inhibits its filament association.
Biochem Biophys Res Commun 477, 820–825.



**Tretyakova I et al. (2016).**

Preparation of quadri-subtype influenza virus-like particles using bovine immunodeficiency virus gag protein.
Virology 487, 163–171.

**Yan X et al. (2016).**

Lack of VEGFR2 signaling causes maldevelopment of the intestinal microvasculature and facilitates necrotizing enterocolitis in neonatal mice.
Am J Physiol Gastrointest Liver Physiol 310, G716–G725.

**Zhou X et al. (2016).**

Cecropin B represses CYP3A29 expression through activation of the TLR2/4-NF- κ B/PXR signaling pathway.
Sci Rep 6, 27876.

**Dong, X et al. (2015).**

BmREEPa is a novel gene that facilitates BmNPV entry into silkworm cells.
PLoS One 10, e0144575.

**Kaiser A et al. (2015).**

A putative non-canonical Ras-like GTPase from *P. falciparum*: chemical properties and characterization of the protein.
PLoS One 10, e0140994.

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

Activation of CXCL-8 transcription by Hepatitis E virus ORF-1 via AP-1.
Mediators Inflamm 2015, 495370.

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

Osteoprotegerin induces apoptosis of osteoclasts and osteoclast precursor cells via the Fas/Fas ligand pathway.
PLoS One 10, e0142519.

**Menden H et al. (2015).**

Lipopolysaccharide (LPS)-mediated angiotensin-2-dependent autocrine angiogenesis is regulated by NADPH oxidase 2 (Nox2) in human pulmonary microvascular endothelial cells.
J Biol Chem 290, 5,449–5,461.



Visit bio-rad.com/getsurebeads for more information.

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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

