

# TGX™ and TGX Stain-Free™ Gels: Publications List



TGX Stain-Free Gels

Bulletin 6507

## 2013

### **De Petrocellis L et al. (2013).**

Non-THC cannabinoids inhibit prostate carcinoma growth in vitro and in vivo: Pro-apoptotic effects and underlying mechanisms. *Br J Pharmacol* 168, 79–102.

---

### **Gallo A et al. (2013).**

Gross cystic disease fluid protein-15(GCDFP-15)/prolactin-inducible protein (PIP) as functional salivary biomarker for primary Sjögren's syndrome. *J Genet Syndr Gene Ther* 4.

---

### **Gan HT et al. (2013).**

Characterization and removal of aggregates formed by nonspecific interaction of IgM monoclonal antibodies with chromatin catabolites during cell culture production. *J Chromatogr A* 1291, 33–40.

---

### **Hegerle N et al. (2013).**

In-vitro and in-vivo analysis of the production of the bordetella type three secretion system effector a in bordetella pertussis, bordetella parapertussis and bordetella bronchiseptica. *Microbes Infect* 15, 399–408.

---

### **Kanzaki H et al. (2013).**

The keap1/nrf2 protein axis plays a role in osteoclast differentiation by regulating intracellular reactive oxygen species signaling. *J Biol Chem* 288, 23009–23020.

---

### **Niemi P et al. (2013).**

Pre-hydrolysis with carbohydrases facilitates the release of protein from brewer's spent grain. *Bioresour Technol* 136, 529–534.

---

### **Parviainen VI et al. (2013).**

Label-free mass spectrometry proteome quantification of human embryonic kidney cells following 24 hours of sialic acid overproduction. *Proteome Sci* 11, 38.

---

### **Rainczuk A et al. (2013).**

The utility of isotope-coded protein labeling for prioritization of proteins found in ovarian cancer patient urine. *J Proteome Res* 12, 4074–4088.

---

**BIO-RAD**

**Roach G et al. (2013).**

Loss of *ascl1a* prevents secretory cell differentiation within the zebrafish intestinal epithelium resulting in a loss of distal intestinal motility. *Dev Biol* 376, 171–186.

---

**Salvachua D et al. (2013).**

Versatile peroxidase as a valuable tool for generating new biomolecules by homogeneous and heterogeneous cross-linking. *Enzyme Microb Technol* 52, 303–311.

---

**Skovgaard PA and Jorgensen H (2013).**

Influence of high temperature and ethanol on thermostable lignocellulolytic enzymes. *J Ind Microbiol Biotechnol* 40, 447–456.

---

**Warmerdam A et al. (2013).**

Characterization of  $\beta$ -galactosidase isoforms from *Bacillus circulans* and their contribution to GOS production. *Appl Biochem Biotechnol* 170, 340–358.

---

**2012**

**Dutka TL et al. (2012).**

Effects of carnosine on contractile apparatus  $Ca^{2+}$  sensitivity and sarcoplasmic reticulum  $Ca^{2+}$  release in human skeletal muscle fibers. *J Appl Physiol* 112, 728–736.

---

**Eckert C et al. (2012).**

Genetic analysis of the Hox hydrogenase in the cyanobacterium *Synechocystis* sp. PCC 6803 reveals subunit roles in association, assembly, maturation, and function. *J Biol Chem* 287, 43502–43515.

---

**Hassiotou F et al. (2012).**

Breastmilk is a novel source of stem cells with multilineage differentiation potential. *Stem Cells* 30, 2164–2174.

---

**Lee J et al. (2012).**

Principles and applications of steric exclusion chromatography. *J Chromatogr A* 1270, 162–170.

---

**Mollica JP et al. (2012).**

S-glutathionylation of troponin i (fast) increases contractile apparatus  $Ca^{2+}$  sensitivity in fast-twitch muscle fibres of rats and humans. *J Physiol* 590, 1443–1463.

---

**Peiris H et al. (2012).**

Increased expression of the glucose-responsive gene, RCAN1, causes hypoinsulinemia,  $\beta$ -cell dysfunction, and diabetes. *Endocrinology* 153, 5212–5221.

---

**Sanata OK et al. (2012).**

Treatments with xylanase at high (90%) and low (40%) water content have different impacts on physicochemical properties of wheat bran. *Food Bioprocess Technol* 10, published online ahead of print October, 2012, accessed September 24, 2013.

---

## TGX Gels

### 2013

**Alconcel SN et al. (2013).**

Synthesis of biotinylated aldehyde polymers for biomolecule conjugation. *Macromol Rapid Commun* 34, 983–989.

---

**Ando H et al. (2013).**

A hypoxia-inducible factor (HIF)-3 $\alpha$  splicing variant, HIF-3 $\alpha$ 4 impairs angiogenesis in hypervascular malignant meningiomas with epigenetically silenced HIF-3 $\alpha$ 4. *Biochem Biophys Res Commun* 433, 139–144.

---

**Apolloni S et al. (2013).**

The NADPH oxidase pathway is dysregulated by the P2X7 receptor in the SOD1-G93A microglia model of amyotrophic lateral sclerosis. *J Immunol* 190, 5187–5195.

---

**Baldini C et al. (2013).**

The P2X receptor-inflammasome complex has a role in modulating the inflammatory response in primary Sjogren's syndrome. *J Intern Med*, published online ahead of print August 1, 2013, accessed September 24, 2013.

---

**Bengtsson AM et al. (2013).**

The cysteinyl leukotriene 2 receptor contributes to all-trans retinoic acid-induced differentiation of colon cancer cells. *BMC Cancer* 13, 336.

---

**Bovell AM and Warncke K (2013).**

The structural model of *Salmonella typhimurium* ethanolamine ammonia-lyase directs a rational approach to the assembly of the functional [(EutB-EutC)<sub>2</sub>]<sub>3</sub> oligomer from isolated subunits. *Biochemistry* 52, 1419–1428.

---

**Brindley MA et al. (2013).**

A stabilized headless measles virus attachment protein stalk efficiently triggers membrane fusion. *J Virol*, published online ahead of print August 21, 2013, accessed September 24, 2013.

---

**Caballero A et al. (2013).**

Differential regulation of parvalbumin and calretinin interneurons in the prefrontal cortex during adolescence. *Brain Struct Funct*, published online ahead of print February 12, 2013, accessed September 24, 2013.

---

**Carbone JW et al. (2013).**

Effects of energy deficit, dietary protein, and feeding on intracellular regulators of skeletal muscle proteolysis. *FASEB J*, published online ahead of print August 21, 2013, accessed September 24, 2013.

---

**Ceglia L et al. (2013).**

Effects of alkali supplementation and vitamin D insufficiency on rat skeletal muscle. *Endocrine*, published online ahead of print May 11, 2013, accessed September 24, 2013.

---

**Chattopadhyay A et al. (2013).**

A novel approach to oral apoA-I mimetic therapy. *J Lipid Res* 54, 995–1010.

---

**Chechenova MB et al. (2013).**

The *Drosophila* Z-disc protein Z(210) is an adult muscle isoform of Zasp52, which is required for normal myofibril organization in indirect flight muscles. *J Biol Chem* 288, 3718–3726.

---

**Chen J et al. (2013).**

The large groove found in the gH/gL structure is an important functional domain for Epstein-Barr virus fusion. *J Virol* 87, 3620–3627.

---

**Chen J et al. (2013).**

5'-UTR mediated translational control of splicing assembly factor RNP-4F expression during development of the *Drosophila* central nervous system. *Gene* 528, 154–162.

---

**di Martino E et al. (2013).**

An integrated genomic, transcriptional and protein investigation of FGFR1 as a putative 4p16.3 deletion target in bladder cancer. *Genes Chromosomes Cancer* 52, 860–871.

---

**Dodd RD et al. (2013).**

NF1 deletion generates multiple subtypes of soft-tissue sarcoma that respond to MEK inhibition. *Mol Cancer Ther* 12, 1906–1917.

---

**Edwards IJ et al. (2013).**

Na<sup>+</sup>/K<sup>+</sup> ATPase  $\alpha$ 1 and  $\alpha$ 3 isoforms are differentially expressed in  $\alpha$ - and  $\gamma$ -motoneurons. *J Neurosci* 33, 9913–9919.

---

**Evers MM et al. (2013).**

Ataxin-3 protein modification as a treatment strategy for spinocerebellar ataxia type 3: Removal of the CAG containing exon. *Neurobiol Dis* 58, 49–56.

---

**Fagan-Solis KD et al. (2013).**

The RhoA pathway mediates MMP-2 and MMP-9-independent invasive behavior in a triple-negative breast cancer cell line. *J Cell Biochem* 114, 1385–1394.

---

**Fischer I et al. (2013).**

Continuous protein purification using functionalized magnetic nanoparticles in aqueous micellar two-phase systems. *J Chromatogr A* 1305, 7–16.

---

**Francin-Allami M et al. (2013).**

Comparative study of wheat low-molecular-weight glutenin and  $\alpha$ -gliadin trafficking in tobacco cells. *Plant Cell Rep* 32, 89–101.

---

**Francis SP et al. (2013).**

A novel role of cytosolic protein synthesis inhibition in aminoglycoside ototoxicity. *J Neurosci* 33, 3079–3093.

---

**Fraser KB et al. (2013).**

LRRK2 secretion in exosomes is regulated by 14-3-3. *Hum Mol Genet*, published online ahead of print August 6, 2013, accessed September 24, 2013.

---

**Fu Y et al (2013).**

An internal standard for protein purification by affinity sorption. *Anal Methods* 5, 1352–1354.

---

**Gee EP et al. (2013).**

SLGISWD sequence in the 10FNIII domain initiates fibronectin fibrillogenesis. *J Biol Chem* 288, 21329–21340.

---

**Hamzic N et al. (2013).**

Immune-induced expression of lipocalin-2 in brain endothelial cells: Relationship with interleukin-6, cyclooxygenase-2 and the febrile response. *J Neuroendocrinol* 25, 271–280.

---

**Harrison LM et al. (2013).**

Effects of the RAS homolog Rhes on Akt/protein kinase B and glycogen synthase kinase 3 phosphorylation in striatum. *Neuroscience* 236, 21–30.

---

**Hasim S et al. (2013).**

Histone biotinylation in *Candida albicans*. *FEMS Yeast Res* 13, 529–539.

---

**Hebert SL et al. (2013).**

The role of pitx2 in maintaining the phenotype of myogenic precursor cells in the extraocular muscles. *PLoS One* 8, e58405.

---

**Heckler EJ et al. (2013).**

Protein disulfide-isomerase interacts with soluble guanylyl cyclase via a redox-based mechanism and modulates its activity. *Biochem J* 452, 161–169.

---

**Hilden K et al. (2013).**

Heterologous expression and structural characterization of two low pH laccases from a biopulping white-rot fungus *Physisporinus rivulosus*. *Appl Microbiol Biotechnol* 97, 1589–1599.

---

**Johnson AC et al. (2013).**

A mutation in the start codon of  $\gamma$ -crystallin D leads to nuclear cataracts in the Dahl SS/Jr-Ctr strain. *Mamm Genome* 24, 95–104.

---

**Kase ET et al. (2013).**

Remodeling of oxidative energy metabolism by galactose improves glucose handling and metabolic switching in human skeletal muscle cells. *PLoS One* 8, e59972.

---

**Kawada I et al. (2013).**

Paxillin mutations affect focal adhesions and lead to altered mitochondrial dynamics: Relevance to lung cancer. *Cancer Biol Ther* 14, 679–691.

---

**Kedees MH et al. (2013).**

Functional activity of murine intestinal mucosal cells is regulated by the glucagon-like peptide-1 receptor. *Peptides* 48C, 36–44.

---

**Khan S et al. (2013).**

MiR-379 regulates cyclin B1 expression and is decreased in breast cancer. *PLoS One* 8, e68753.

---

**Kinney MA et al. (2013).**

Temporal modulation of  $\beta$ -catenin signaling by multicellular aggregation kinetics impacts embryonic stem cell cardiomyogenesis. *Stem Cells Dev* 22, 2665–2677.

---

**Kirby ED et al. (2013).**

Acute stress enhances adult rat hippocampal neurogenesis and activation of newborn neurons via secreted astrocytic FGF2. *Elife* 2, e00362.

---

**Kjaer TR et al. (2013).**

M-ficolin binds selectively to the capsular polysaccharides of *Streptococcus pneumoniae* serotypes 19B and 19C and of a *Streptococcus mitis* strain. *Infect Immun* 81, 452–459.

---

**Komiyama H et al. (2013).**

Cyclical mechanical stretch enhances degranulation and IL-4 secretion in RBL-2H3 mast cells. *Cell Biochem Funct*, published online ahead of print April 15, 2013, accessed September 24, 2013.

---

**Kumar S et al. (2013).**

Conformational changes of recombinant Ca(2+) -ATPase studied by reaction-induced infrared difference spectroscopy. *FEBS J*.

---

**Kuramitsu Y et al. (2013).**

Up-regulation of ddx39 in human malignant pleural mesothelioma cell lines compared to normal pleural mesothelial cells. *Anticancer Res* 33, 2557–2560.

---

**Langevin C et al. (2013).**

Zebrafish ISG15 exerts a strong antiviral activity against RNA and DNA viruses and regulates the interferon response. *J Virol* 87, 10025–10036.

---

**Lauer ME et al. (2013).**

Irreversible heavy chain transfer to hyaluronan oligosaccharides by tumor necrosis factor-stimulated gene-6. *J Biol Chem* 288, 205–214.

---

**Lauer ME et al. (2013).**

Tumor necrosis factor-stimulated gene-6 (TSG-6) amplifies hyaluronan synthesis by airway smooth muscle cells. *J Biol Chem* 288, 423–431.

---

**Leclere R et al. (2013).**

Effect of estrogens on base excision repair in brain and liver mitochondria of aged female rats. *Biogerontology* 14, 383–394.

---

**Li L et al. (2013).**

Parvoviral left-end hairpin ears are essential during infection for establishing a functional intranuclear transcription template and for efficient progeny genome encapsidation. *J Virol* 87, 10501–10514.

---

**Liu B et al. (2013).**

Astrocyte elevated gene-1 regulates osteosarcoma cell invasion and chemoresistance via endothelin-1/endothelin A receptor signaling. *Oncol Lett* 5, 505–510.

---

**Long AR et al. (2013).**

A detergent-free strategy for the reconstitution of active enzyme complexes from native biological membranes into nanoscale discs. *BMC Biotechnol* 13, 41.

---

**Lu J et al. (2013).**

Bradykinin B2 receptor contributes to the exaggerated muscle mechanoreflex in rats with femoral artery occlusion. *Am J Physiol Heart Circ Physiol* 304, H1166–1174.

---

**Luo H et al. (2013).**

Resveratrol induces premature senescence in lung cancer cells via ROS-mediated DNA damage. *PLoS One* 8, e60065.

---

**Maire V et al. (2013).**

TTK/hMPS1 is an attractive therapeutic target for triple-negative breast cancer. *PLoS One* 8, e63712.

---

**Marshall AC et al. (2013).**

Fetal betamethasone exposure attenuates angiotensin-(1-7)-Mas receptor expression in the dorsal medulla of adult sheep. *Peptides* 44, 25–31.

---

**Martinez-Pastor M et al. (2013).**

Negative feedback regulation of the yeast CTH1 and CTH2 mRNA binding proteins is required for adaptation to iron deficiency and iron supplementation. *Mol Cell Biol* 33, 2178–2187.

---

**Martorell P et al. (2013).**

A cocoa peptide protects *Caenorhabditis elegans* from oxidative stress and  $\beta$ -amyloid peptide toxicity. *PLoS One* 8, e63283.

---

**Matsumoto NM et al. (2013).**

Synthesis of nanogel–protein conjugates. *Polym Chem* 84, 2464–2469.

---

**Matsumoto NM et al. (2013).**

Smart vaults: Thermally-responsive protein nanocapsules. *ACS Nano* 7, 867–874.

---

**Meng QX et al. (2013).**

Association of serum angiopoietin-like protein 2 and epinephrine levels in metabolically healthy but obese individuals: In vitro and in vivo evidence. *Exp Ther Med* 5, 1631–1636.

---

**Miyake M et al. (2013).**

Expression of CXCL1 in human endothelial cells induces angiogenesis through the CXCR2 receptor and the ERK1/2 and EGF pathways. *Lab Invest* 93, 768–778.

---

**Moore CJ et al. (2013).**

Malondialdehyde and 4-hydroxynonenal adducts are not formed on cardiac ryanodine receptor (RyR2) and sarco(endo)plasmic reticulum Ca<sup>2+</sup>-ATPase (SERCA2) in diabetes. *Mol Cell Biochem* 376, 121–135.

---

**Moy GA and McNay EC (2013).**

Caffeine prevents weight gain and cognitive impairment caused by a high-fat diet while elevating hippocampal BDNF. *Physiol Behav* 109, 69–74.

---

**Nair S et al. (2013).**

The chromosomal passenger protein birc5b organizes microfilaments and germ plasm in the zebrafish embryo. *PLoS Genet* 9, e1003448.

---

**Neman J et al. (2013).**

Co-evolution of breast-to-brain metastasis and neural progenitor cells. *Clin Exp Metastasis*, published online ahead of print February 28, 2013, accessed September 24, 2013.

---

**Nikam A et al. (2013).**

Transition between acute and chronic hepatotoxicity in mice is associated with impaired energy metabolism and induction of mitochondrial heme oxygenase-1. *PLoS One* 8, e66094.

---

**Nohata N et al. (2013).**

Tumour-suppressive microRNA-874 contributes to cell proliferation through targeting of histone deacetylase 1 in head and neck squamous cell carcinoma. *Br J Cancer* 108, 1648–1658.

---

**Nomura Y et al. (2013).**

Reduction of E-cadherin by human defensin-5 in esophageal squamous cells. *Biochem Biophys Res Commun* 439, 71–77.

---

**O'Dell LE et al. (2013).**

Enhanced nicotine self-administration and suppressed dopaminergic systems in a rat model of diabetes. *Addict Biol*, epub ahead of print.

---

**O'Farrell F et al. (2013).**

Two-tiered control of epithelial growth and autophagy by the insulin receptor and the ret-like receptor, stitcher. *PLoS Biol* 11, e1001612.

---

**Ohta K et al. (2013).**

Depletion of JARID1B induces cellular senescence in human colorectal cancer. *Int J Oncol* 42, 1212–1218.

---



**Overcash RF et al. (2013).**

Androgen signaling promotes translation of TMEFF2 in prostate cancer cells via phosphorylation of the alpha subunit of the translation initiation factor 2. *PLoS One* 8, e55257.

---

**Oyola-Robles D et al. (2013).**

Identification of novel protein domains required for the expression of an active dehydratase fragment from a polyunsaturated fatty acid synthase. *Protein Sci* 22, 954–963.

---

**Pancani T et al. (2013).**

Effect of high-fat diet on metabolic indices, cognition, and neuronal physiology in aging G344 rats. *Neurobiol Aging* 34, 1977–1987.

---

**Paz EA et al. (2013).**

Polyamines are on cometabolites that regulate the LIN28/let-7 pathway in colorectal cancer cells. *Mol Carcinog*, epub ahead of print.

---

**Qi LM and Wade J (2013).**

Sexually dimorphic and developmentally regulated expression of tubulin-specific chaperone protein A in the LMAN of zebra finches. *Neuroscience* 247, 182–190.

---

**Rahn JJ et al. (2013).**

Opa1 is required for proper mitochondrial metabolism in early development. *PLoS One* 8, e59218.

---

**Restaino OF et al. (2013).**

High cell density cultivation of a recombinant *E. coli* strain expressing a key enzyme in bioengineered heparin production. *Appl Microbiol Biotechnol* 97, 3893–3900.

---

**Rogers SC et al. (2013).**

Exposure to high or low glucose levels accelerates the appearance of markers of endothelial cell senescence and induces dysregulation of nitric oxide synthase. *J Gerontol A Biol Sci Med Sci*, published online ahead of print April 12, 2013, accessed September 24, 2013.

---

**Rustagi A et al. (2013).**

Two new monoclonal antibodies for biochemical and flow cytometric analyses of human interferon regulatory factor-3 activation, turnover, and depletion. *Methods* 59, 225–232.

---

**Sakurai T et al. (2013).**

Cardiomyocyte FGF signaling is required for Cx43 phosphorylation and cardiac gap junction maintenance. *Exp Cell Res* 319, 2152–2165.

---

**Scott DW et al. (2013).**

Heterogenic endothelial responses to inflammation: Role for differential N-glycosylation and vascular bed of origin. *J Am Heart Assoc* 2, e000263.

---

**Scott DW et al. (2013).**

Identification of a high-mannose ICAM-1 glycoform: Effects of ICAM-1 hypoglycosylation on monocyte adhesion and outside in signaling. *Am J Physiol Cell Physiol* 305, C228–237.

---

**Shavlakadze T et al. (2013).**

Short-term feed deprivation rapidly induces the protein degradation pathway in skeletal muscles of young mice. *J Nutr* 143, 403–409.

---

**Shavlakadze T et al. (2013).**

Impact of fasting on the rhythmic expression of myogenic and metabolic factors in skeletal muscle of adult mice. *Am J Physiol Cell Physiol* 305, C26–35.

---

**Swaidani S et al. (2013).**

TSG-6 protein is crucial for the development of pulmonary hyaluronan deposition, eosinophilia, and airway hyperresponsiveness in a murine model of asthma. *J Biol Chem* 288, 412–422.

---

**Tanca A et al. (2013).**

Comparison of detergent-based sample preparation workflows for LTQ-Orbitrap analysis of the *Escherichia coli* proteome. *Proteomics* 13, 2597–2607

---

**Tang AT et al. (2013).**

ROCK1 feedback regulation of the upstream small GTPase RhoA. *Cell Signal* 24, 1375–1380.

---

**Tang-Feldman YJ et al. (2013).**

Murine cytomegalovirus (MCMV) infection upregulates P38 MAP kinase in aortas of Apo E KO mice: A molecular mechanism for MCMV-induced acceleration of atherosclerosis. *J Cardiovasc Transl Res* 6, 54–64.

---

**Troxell B et al. (2013).**

Manganese and zinc regulate virulence determinants in *Borrelia burgdorferi*. *Infect Immun* 81, 2743–2752.

---

**Walker EJ et al. (2013).**

Rhinovirus 3C protease facilitates specific nucleoporin cleavage and mislocalisation of nuclear proteins in infected host cells. *PLoS One* 8, e71316.

---

**Weekley CM et al. (2013).**

Selenium metabolism in cancer cells: The combined application of XAS and XFM techniques to the problem of selenium speciation in biological systems. *Nutrients* 5, 1734–1756.

---

**Wijayanti HB et al. (2013).**

Reduction of aggregation of  $\beta$ -lactoglobulin during heating by dihydrolipoic acid. *J Dairy Res* 1–7.

---

**Yamamoto N et al. (2013).**

Tumor suppressive microRNA-218 inhibits cancer cell migration and invasion by targeting focal adhesion pathways in cervical squamous cell carcinoma. *Int J Oncol* 42, 1523–1532.

---

**Yap MN and Bernstein HD (2013).**

Mutations in the *Escherichia coli* ribosomal protein L22 selectively suppress the expression of a secreted bacterial virulence factor. *J Bacteriol* 195, 2991–2999.

---

**Yoshioka Y et al. (2013).**

Comparative marker analysis of extracellular vesicles in different human cancer types. *J Extracell Vesicles* 2, ev.v2i0.20424.

---

**Yousef H et al. (2013).**

hESC-secreted proteins can be enriched for multiple regenerative therapies by heparin-binding. *Aging (Albany NY)* 5, 357–372.

---

**Zhang L et al. (2013).**

Deletion of *Synechocystis* sp. PCC 6803 leader peptidase LepB1 affects photosynthetic complexes and respiration. *Mol Cell Proteomics* 12, 1192–1203.

---

**Zhou Y et al. (2013).**

TWIST interacts with endothelin-1/endothelin A receptor signaling in osteosarcoma cell survival against cisplatin. *Oncol Lett* 5, 857–861.

---

## 2012

**Arumugam A et al. (2012).**

Mammary cancer promotion by ovarian hormones involves IGFR/AKT/mTOR signaling. *Steroids* 77, 791–797.

---

**Bakkebo M et al. (2012).**

SARA is dispensable for functional TGF- $\beta$  signaling. *FEBS Lett* 586, 3367–3372.

---

**Coltharp C et al. (2012).**

Accurate construction of photoactivated localization microscopy (PALM) images for quantitative measurements. *PLoS One* 7, e51725.

---

**Cully TR et al. (2012).**

Changes in plasma membrane Ca-ATPase and stromal interacting molecule 1 expression levels for Ca(2+) signaling in dystrophic mdx mouse muscle. *Am J Physiol Cell Physiol* 303, C567-576.

---

**Cunningham MD et al. (2012).**

Wapl antagonizes cohesin binding and promotes polycomb-group silencing in *Drosophila*. *Development* 139, 4172–4179.

---

**Degn SE et al. (2012).**

Mannan-binding lectin-associated serine protease (MASP)-1 is crucial for lectin pathway activation in human serum, whereas neither MASP-1 nor MASP-3 is required for alternative pathway function. *J Immunol* 189, 3957–3969.

---

**De Leo A et al. (2012).**

Resveratrol inhibits Epstein Barr virus lytic cycle in Burkitt's lymphoma cells by affecting multiple molecular targets. *Antiviral Res* 96, 196–202.

---

**Delobel J et al. (2012).**

Subcellular fractionation of stored red blood cells reveals a compartment-based protein carbonylation evolution. *J Proteomics* 76 Spec No., 181–193.

---

**D'Orso I et al. (2012).**

Transition step during assembly of HIV Tat:P-TEFb transcription complexes and transfer to TAR RNA. *Mol Cell Biol* 32, 4780–4793

---

**Fischer I and Franzreb M (2012).**

Nanoparticle mediated protein separation in aqueous micellar two-phase systems. *Solvent Extraction and Ion Exchange* 30, 1–16.

---

**Franzetti E et al. (2012).**

Autophagy precedes apoptosis during the remodeling of silkworm larval midgut. *Apoptosis* 17, 305–324.

---

**Gabery S et al. (2012).**

Characterization of a rat model of Huntington's disease based on targeted expression of mutant huntingtin in the forebrain using adeno-associated viral vectors. *Eur J Neurosci* 36, 2789–2800.

---

**Gao M et al. (2012).**

The role of forkhead box q1 transcription factor in ovarian epithelial carcinomas. *Int J Mol Sci* 13, 13881–13893.

---

**Gestl EE and Anne Bottger S (2012).**

Cytoplasmic sequestration of the tumor suppressor p53 by a heat shock protein 70 family member, mortalin, in human colorectal adenocarcinoma cell lines. *Biochem Biophys Res Commun* 423, 411–416.

---

**Gibson MS et al. (2012).**

Identification, cloning, and functional characterization of the IL-1 receptor antagonist in the chicken reveal important differences between the chicken and mammals. *J Immunol* 189, 539–550.

---

**Gilbert NM et al. (2012).**

A glycosylphosphatidylinositol anchor is required for membrane localization but dispensable for cell wall association of chitin deacetylase 2 in *Cryptococcus neoformans*. *MBio* 3, e00007-12

---

**Gnosa S et al. (2012).**

Expression of AEG-1 mRNA and protein in colorectal cancer patients and colon cancer cell lines. *J Transl Med* 10, 109.

---

**Hardison SE et al. (2012).**

Protective immunity against pulmonary cryptococcosis is associated with STAT1-mediated classical macrophage activation. *J Immunol* 189, 4060–4068.

---

**Hare JM et al. (2012).**

The *Acinetobacter* regulatory UmuDAb protein cleaves in response to DNA damage with chimeric LexA/UmuD characteristics. *FEMS Microbiol Lett* 334, 57–65.

---

**Hess K et al. (2012).**

Effects of MASP-1 of the complement system on activation of coagulation factors and plasma clot formation. *PLoS One* 7, e35690.

---

**Hessvik NP et al. (2012).**

Profiling of microRNAs in exosomes released from PC-3 prostate cancer cells. *Biochim Biophys Acta* 1819, 1154–1163.

---

**Hughes, JB et al. (2012).**

Pertuzumab increases 17-AAG-induced degradation of ErbB2, and this effect is further increased by combining pertuzumab with trastuzumab. *Pharmaceuticals* 5, 674–689.

---

**Ikenaga N et al. (2012).**

Pancreatic cancer cells enhance the ability of collagen internalization during epithelial-mesenchymal transition. *PLoS One* 7, e40434.

---

**Isozaki Y et al. (2012).**

Identification of novel molecular targets regulated by tumor suppressive miR-375 induced by histone acetylation in esophageal squamous cell carcinoma. *Int J Oncol* 41, 985–994.

---

**Jarome TJ et al. (2012).**

The timing of multiple retrieval events can alter GluR1 phosphorylation and the requirement for protein synthesis in fear memory reconsolidation. *Learn Mem* 19, 300–306.

---

**Kim S et al. (2012).**

Efficacy of phosphatidylinositol-3 kinase inhibitors in a primary mouse model of undifferentiated pleomorphic sarcoma. *Sarcoma* 2012, 680708.

---

**Kinoshita T et al. (2012).**

Tumor suppressive microRNA-375 regulates lactate dehydrogenase B in maxillary sinus squamous cell carcinoma. *Int J Oncol* 40, 185–193.

---

**Kinoshita T et al. (2012).**

Tumor suppressive microRNA-133a regulates novel targets: Moesin contributes to cancer cell proliferation and invasion in head and neck squamous cell carcinoma. *Biochem Biophys Res Commun* 418, 378–383.

---

**Kinoshita T et al. (2012).**

Actin-related protein 2/3 complex subunit 5 (ARPC5) contributes to cell migration and invasion and is directly regulated by tumor-suppressive microRNA-133a in head and neck squamous cell carcinoma. *Int J Oncol* 40, 1770–1778.

---

**Kobayashi R et al. (2012).**

Trichloroethylene enhances TCR-CD3-induced proliferation of CD8(+) rather than CD4(+) T cells. *J Toxicol Sci* 37, 381–387.

---

**Krammer F et al. (2012).**

A carboxy-terminal trimerization domain stabilizes conformational epitopes on the stalk domain of soluble recombinant hemagglutinin substrates. *PLoS One* 7, e43603.

---

**Kumaresan A et al. (2012).**

Oviductal fluid modulates the dynamics of tyrosine phosphorylation in cryopreserved boar spermatozoa during capacitation. *Mol Reprod Dev* 79, 525–540.

---

**Kumaresan A et al. (2012).**

Quantification of kinetic changes in protein tyrosine phosphorylation and cytosolic Ca<sup>2+</sup> concentration in boar spermatozoa during cryopreservation. *Reprod Fertil Dev* 24, 531–542.

---

**Lappas M (2012).**

Anti-inflammatory properties of sirtuin 6 in human umbilical vein endothelial cells. *Mediators Inflamm* 2012, 597514.

---

**Lee JJ et al. (2012).**

Seroproteomics reveals cross-reactivity of anti-enterovirus 71 antibodies with cytoplasmic actin. *International Journal of Virology and Molecular Biology* 2, 12–17

---

**Li J and Murtaugh MP (2012).**

Dissociation of porcine reproductive and respiratory syndrome virus neutralization from antibodies specific to major envelope protein surface epitopes. *Virology* 433, 367–376

---

**Liu X et al. (2012).**

Heterogeneous expression of the core circadian clock proteins among neuronal cell types in mouse retina. *PLoS One* 7, e50602.

---

**McCaffrey G et al. (2012).**

P-glycoprotein trafficking at the blood-brain barrier altered by peripheral inflammatory hyperalgesia. *J Neurochem* 122, 962–975.

---

**McGreal RS et al. (2012).**

$\alpha\beta$ -crystallin/sHSP protects cytochrome c and mitochondrial function against oxidative stress in lens and retinal cells. *Biochim Biophys Acta* 1820, 921–930.

---

**Miyoshi J et al. (2012).**

5-aminosalicylic acid mediates expression of cyclooxygenase-2 and 15-hydroxyprostaglandin dehydrogenase to suppress colorectal tumorigenesis. *Anticancer Res* 32, 1193–1202.

---

**Monteiro KM et al. (2012).**

*Echinococcus granulosus* antigen B structure: Subunit composition and oligomeric states. *PLoS Negl Trop Dis* 6, e1551.

---

**Musha A et al. (2012).**

Synergistic effect of heat shock protein 90 inhibitor, 17-allylamino-17-demethoxygeldanamycin and X-rays, but not carbon-ion beams, on lethality in human oral squamous cell carcinoma cells. *J Radiat Res* 53, 545–550.

---

**Muttray A et al. (2012).**

Haemocytic leukemia in Prince Edward Island (PEI) soft shell clam (*Mya arenaria*): Spatial distribution in agriculturally impacted estuaries. *Sci Total Environ* 424, 130–142.

---

**Natabotu SK et al. (2012).**

Control of myosin motor activity by the reversible alteration of protein structure for application in a bionanodevice. *Adv Sci Lett* 16, 213–221.

---

**Patankar JV et al. (2012).**

Intestinal GATA4 deficiency protects from diet-induced hepatic steatosis. *J Hepatol* 57, 1061–1068.

---

**Rabadi MM et al. (2012).**

Interaction between uric acid and HMGB1 translocation and release from endothelial cells. *Am J Physiol Renal Physiol* 302, F730–741.

---

**Rabadi MM et al. (2012).**

HMGB1 in renal ischemic injury. *Am J Physiol Renal Physiol* 303, F873–885.

---

**Ratchford SM et al. (2012).**

Proteins regulating cap-dependent translation are downregulated during total knee arthroplasty. *Am J Physiol Regul Integr Comp Physiol* 302, R702–711.

---

**Rommelfanger DM et al. (2012).**

Systemic combination virotherapy for melanoma with tumor antigen-expressing vesicular stomatitis virus and adoptive T-cell transfer. *Cancer Res* 72, 4753–4764.

---

**Simmons S et al. (2012).**

Biphenotypic B-lymphoid/myeloid cells expressing low levels of Pax5: Potential targets of BAL development. *Blood* 120, 3688–3698.

---

**Tanca A et al. (2012).**

Comparability of differential proteomics data generated from paired archival fresh-frozen and formalin-fixed samples by GeLC-MS/MS and spectral counting. *J Proteomics* 77, 561–576.

---

**Tanimoto Y et al. (2012).**

Prevention of premature fusion of calvarial suture in GLI-Kruppel family member 3 (Gli3)-deficient mice by removing one allele of Runt-related transcription factor 2 (Runx2). *J Biol Chem* 287, 21429–21438.

---

**Tenayuca JM and Nazarian A (2012).**

Hydrocodone and morphine possess similar rewarding effects and reduce ERK and CREB phosphorylation in the nucleus accumbens. *Synapse* 66, 918–922.

---

**Tetreau G et al. (2012).**

Larval midgut modifications associated with Bti resistance in the yellow fever mosquito using proteomic and transcriptomic approaches. *BMC Genomics* 13, 248.

---

**Vishwanatha KS et al. (2012).**

Structural organization of the nine spectrin repeats of Kalirin. *Biochemistry* 51, 5663–5673.

---

**Wan YW et al. (2012).**

A smoking-associated 7-gene signature for lung cancer diagnosis and prognosis. *Int J Oncol* 41, 1387–1396.

---

**Watanabe J et al. (2012).**

Proteomic profiling following immunoaffinity capture of high-density lipoprotein: Association of acute-phase proteins and complement factors with proinflammatory high-density lipoprotein in rheumatoid arthritis. *Arthritis Rheum* 64, 1828–1837.

---

**Weber E et al. (2012).**

GFP facilitates native purification of recombinant perlucin derivatives and delays the precipitation of calcium carbonate. *PLoS One* 7, e46653.

---

**Whitcomb SJ et al. (2012).**

Histone monoubiquitylation position determines specificity and direction of enzymatic cross-talk with histone methyltransferases Dot1L and PRC2. *J Biol Chem* 287, 23718–23725.

---

**Wolschendorf F et al. (2012).**

Kinase control prevents HIV-1 reactivation in spite of high levels of induced NF- $\kappa$ B activity. *J Virol* 86, 4548–4558.

---

**Xing J et al. (2012).**

Angiotensin-(1-7) inhibits neuronal activity of dorsolateral periaqueductal gray via a nitric oxide pathway. *Neurosci Lett* 522, 156–161.

---

**Zou S et al. (2012).**

Notch signaling in descending thoracic aortic aneurysm and dissection. *PLoS One* 7, e52833.

---

## 2011

**Ball KK et al. (2011).**

Reduced gap junctional communication among astrocytes in experimental diabetes: Contributions of altered connexin protein levels and oxidative-nitrosative modifications. *J Neurosci Res* 89, 2052–2067.

---

**Chu VC et al. (2011).**

Persistent expression of hepatitis C virus non-structural proteins leads to increased autophagy and mitochondrial injury in human hepatoma cells. *PLoS One* 6, e28551.

---

**Kallberg K et al. (2011).**

Application of a pH responsive multimodal hydrophobic interaction chromatography medium for the analysis of glycosylated proteins. *J Chromatogr A* 1218, 678–683.

---

**Kumaresan A et al. (2011).**

Cryopreservation-induced alterations in protein tyrosine phosphorylation of spermatozoa from different portions of the boar ejaculate. *Cryobiology* 63, 137–144.

---



**Monnier PP et al. (2011).**

Involvement of caspase-6 and caspase-8 in neuronal apoptosis and the regenerative failure of injured retinal ganglion cells. *J Neurosci* 31, 10494–10505.

---

**Ngo S and Guo Z (2011).**

Key residues for the oligomerization of A $\beta$ 42 protein in Alzheimer's disease. *Biochem Biophys Res Commun* 414, 512–516.

---

**Ruiz Z et al. (2011).**

Recruitment of DNA replication and damage response proteins to viral replication centers during infection with NS2 mutants of Minute Virus of Mice (MVM). *Virology* 410, 375–384.

---

**Snow JB et al. (2011).**

Role for PKC $\beta$  in enhanced endothelin-1-induced pulmonary vasoconstrictor reactivity following intermittent hypoxia. *Am J Physiol Lung Cell Mol Physiol* 301, L745–754.

---

**Sproul A et al. (2011).**

N-methyl-D-aspartate receptor subunit NR3a expression and function in principal cells of the collecting duct. *Am J Physiol Renal Physiol* 301, F44–54.

---

**Stone DK et al. (2011).**

A model of nitric oxide induced  $\alpha$ -synuclein misfolding in Parkinson's disease. *Neurosci Lett* 523, 167–173.

---

**Yamaguchi K (2011).**

Preparation and proteomic analysis of chloroplast ribosomes. *Methods Mol Biol* 775, 241–264.

---

**Zhang D et al. (2011).**

Establishment of pemetrexed-resistant non-small cell lung cancer cell lines. *Cancer Lett* 309, 228–235.

---

## 2010

**Ambartsumyan G et al. (2010).**

Centromere protein a dynamics in human pluripotent stem cell self-renewal, differentiation and DNA damage. *Hum Mol Genet* 19, 3970–3982.

---



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 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** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000