

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 β -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, β -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 α splicing variant, HIF-3 α 4 impairs angiogenesis in hypervascular malignant meningiomas with epigenetically silenced HIF-3 α 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)₂]₃ 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⁺/K⁺ ATPase α 1 and α 3 isoforms are differentially expressed in α - and γ -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 α -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 γ -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 β -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 β -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 Ca²⁺-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 β -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- β 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 Ca²⁺ 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- κ 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 β 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 β 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 α -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
Laboratories, Inc.**

Life Science
Group

Web site 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