

Molecular Imager<sup>®</sup> VersaDoc<sup>™</sup> System Citations

- Abad MA et al. (2006). Neuronal Pentraxin 1 Contributes to the Neuronal Damage Evoked by Amyloid- $\beta$  and Is Overexpressed in Dystrophic Neurites in Alzheimer's Brain. *J Neurosci* 26, 12735-12747.
- Abram ME and Parniak MA (2005). Virion Instability of Human Immunodeficiency Virus Type 1 Reverse Transcriptase (RT) Mutated in the Protease Cleavage Site between RT p51 and the RT RNase H Domain. *J Virol* 79, 11952-11961.
- Abt M and Rivers DB (2007). Characterization of phenoloxidase activity in venom from the ectoparasitoid *Nasonia vitripennis* (Walker) (Hymenoptera: Pteromalidae). *J Invertebr Pathol* 94, 108-118.
- Ackerman WE, IV et al. (2007). Association of PAT Proteins with Lipid Storage Droplets in Term Fetal Membranes. *Placenta* 28, 465-476.
- Ackerman WE, IV et al. (2004). Epidermal Growth Factor and Interleukin-1 $\beta$  Utilize Divergent Signaling Pathways to Synergistically Upregulate Cyclooxygenase-2 Gene Expression in Human Amnion-Derived WISH Cells. *Biol Reprod* 71, 2079-2086.
- Ackerman WE, IV et al. (2005). Despite Transcriptional and Functional Coordination, Cyclooxygenase-2 and Microsomal Prostaglandin E Synthase-1 Largely Reside in Distinct Lipid Microdomains in WISH Epithelial Cells. *J Histochem Cytochem* 53, 1391-1401.
- Ackerman WE, IV et al. (2005). Modulation of Cytokine-Induced Cyclooxygenase 2 Expression by PPAR $\gamma$  Ligands Through NF $\kappa$ B Signal Disruption in Human WISH and Amnion Cells. *Biol Reprod* 73, 527-535.
- Agler HL et al. (2003). Custom Distinctions in the Interaction of G-protein  $\beta$  Subunits with N-type (CaV2.2) Versus P/Q-type (CaV2.1) Calcium Channels. *J Gen Physiol* 121, 495-510.
- Agostinho P (2007). Role of cyclin-dependent kinase 5 in the neurodegenerative process triggered by amyloid-beta and Prion peptides: Implications for Alzheimer's disease and Prion-related encephalopathies. *Cell Mol Neurobiol* 27, 943-957.
- Aguilar HC et al. (2006). N-Glycans on Nipah Virus Fusion Protein Protect against Neutralization but Reduce Membrane Fusion and Viral Entry. *J Virol* 80, 4878-4889.
- Aguilar HC et al. (2007). Polybasic KKR Motif in the Cytoplasmic Tail of Nipah Virus Fusion Protein Modulates Membrane Fusion by Inside-Out Signaling. *J Virol* 81, 4520-4532.
- Aguirre JI et al. (2007). A Novel Ligand-independent Function of the Estrogen Receptor Is Essential for Osteocyte and Osteoblast Mechanotransduction. *J Biol Chem* 282, 25501-25508.
- Ahmed R and Duncan RF (2004). Translational Regulation of Hsp90 mRNA: aug-proximal 5'-untranslated region elements essential for preferential heat shock translation. *J Biol Chem* 279, 49919-49930.
- Alb JG, Jr. et al. (2007). The pathologies associated with functional titration of phosphatidylinositol transfer protein  $\alpha$  activity in mice. *J Lipid Res* 48, 1857-1872.
- Alvaro AR et al. (2007). NPY in rat retina is present in neurons, in endothelial cells and also in microglial and Muller cells. *Neurochem Int* 50, 757-763.
- Amici C et al. (2006). Herpes Simplex Virus Disrupts NF- $\kappa$ B Regulation by Blocking Its Recruitment to the I $\kappa$ B $\alpha$  Promoter and Directing the Factor on Viral Genes. *J Biol Chem* 281, 7110-7117.
- Amoruso A et al. (2007). Quantification of PPAR- $\gamma$  protein in monocyte/macrophages from healthy smokers and non-smokers: A possible direct effect of nicotine. *Life Sci* 81, 906-915.
- Ann D (2007). Role for KAP1 serine 824 phosphorylation and sumoylation/desumoylation switch in regulating KAP1-mediated transcriptional repression. *J Biol Chem* 282, 36177-36189.
- Auerbach M and Liedtke CM (2007). Role of the scaffold protein RACK1 in apical expression of CFTR. *Am J Physiol Cell Physiol* 293, C294-C304.
- Ayala JE et al. (2007). Chronic Treatment With Sildenafil Improves Energy Balance and Insulin Action in High Fat-Fed Conscious Mice. *Diabetes* 56, 1025-1033.
- Bachmann PS et al. (2005). Dexamethasone resistance in B-cell precursor childhood acute lymphoblastic leukemia occurs downstream of ligand-induced nuclear translocation of the glucocorticoid receptor. *Blood* 105, 2519-2526.
- Bailey LO et al. (2005). The quantification of cellular viability and inflammatory response to stainless steel alloys. *Biomaterials* 26, 5296-5302.
- Bamberger ME et al. (2003). A Cell Surface Receptor Complex for Fibrillar  $\beta$ -Amyloid Mediates Microglial Activation. *J Neurosci* 23, 2665-2674.
- Bannai H et al. (2003). Cloning and Expression of Mitochondrial Heat Shock Protein 70 of *Trypanosoma congolense* and Potential Use as a Diagnostic Antigen. *Clin Diagn Lab Immunol* 10, 926-933.
- Bardales JR et al. (2007). CK2-mediated phosphorylation of a type II regulatory subunit of cAMP-dependent protein kinase from the mollusk *Mytilus galloprovincialis*. *Arch Biochem Biophys* 461, 130-137.
- Barlow PG et al. (2006). The human cationic host defense peptide LL-37 mediates contrasting effects on apoptotic pathways in different primary cells of the innate immune system. *J Leukoc Biol* 80, 509-520.
- Beatty J (2007). The photosystem of *Rhodobacter sphaeroides* assembles with zincbacteriochlorophyll in a bchD (Magnesium chelatase) mutant. *Biochemistry (Mosc)* 46, 12491-12500.
- Beaulieu N et al. (2007). Regulation of RasGRP1 by B Cell Antigen Receptor Requires Cooperativity between Three Domains Controlling Translocation to the Plasma Membrane. *Mol Biol Cell* 18, 3156-3168.
- Belhocine K et al. (2007). Trans-splicing of the LI.LtrB group II intron in *Lactococcus lactis*. *Nucleic Acids Res* 35, 2257-2268.
- Ben Yehuda S et al. (2005). Defining a Centromere-like Element in *Bacillus subtilis* by Identifying the Binding Sites for the Chromosome-Anchoring Protein RacA. *Mol Cell* 17, 773-782.
- Bengtsson L and Wilson KL (2006). Barrier-to-Autointegration Factor Phosphorylation on Ser-4 Regulates Emerin Binding to Lamin A In Vitro and Emerin Localization In Vivo. *Mol Biol Cell* 17, 1154-1163.
- Bennati AM et al. (2006). Sterol dependent regulation of human TM7SF2 gene expression: Role of the encoded 3[beta]-hydroxysterol  $\delta$ 14-reductase in human cholesterol biosynthesis. *Biochimica Biophys Acta (BBA) - Molecular and Cell Biology of Lipids* 1761, 677-685.
- Bergh JJ et al. (2005). Integrin  $\alpha$ V $\beta$ 3 Contains a Cell Surface Receptor Site for Thyroid Hormone that Is Linked to Activation of Mitogen-Activated Protein Kinase and Induction of Angiogenesis. *Endocrinology* 146, 2864-2871.
- Bhadriraju K et al. (2007). Activation of ROCK by RhoA is regulated by cell adhesion, shape, and cytoskeletal tension. *Exp Cell Res* 313, 3616-3623.
- Binek M (2007). Physiological properties and genomic diversity of *Corynebacterium pseudotuberculosis* strain isolates from goats in Poland. *Med Weter* 63, 1474-1477.
- Blaesse P et al. (2006). Oligomerization of KCC2 Correlates with Development of Inhibitory Neurotransmission. *J Neurosci* 26, 10407-10419.

- Blossom SJ and Gilbert KM (2006). Exposure to a Metabolite of the Environmental Toxicant, Trichloroethylene, Attenuates CD4+ T Cell Activation-Induced Cell Death by Metalloproteinase-Dependent FasL Shedding. *Toxicol Sci* 92, 103-114.
- Blossom SJ et al. (2007). Chronic Exposure to a Trichloroethylene Metabolite in Autoimmune-Prone MRL+/+ Mice Promotes Immune Modulation and Alopecia. *Toxicol Sci* 95, 401-411.
- Bocca C et al. (2007). Conjugated linoleic acid inhibits Caco-2 cell growth via ERK-MAPK signaling pathway. *J Nutr Biochem* 18, 332-340.
- Boero S et al. (2007). Modulation of human lung fibroblast functions by ciclosporin: Evidence for its conversion into the active metabolite desisobutyl-ciclosporin. *Immunol Lett* 112, 39-46.
- Bolognesi M et al. (2005). Haeme oxygenase mediates hyporeactivity to phenylephrine in the mesenteric vessels of cirrhotic rats with ascites. *Gut* 54, 1630-1636.
- Bolognesi M et al. (2007). Carbon Monoxide-Mediated Activation of Large-Conductance Calcium-Activated Potassium Channels Contributes to Mesenteric Vasodilatation in Cirrhotic Rats. *J Pharmacol Exp Ther* 321, 187-194.
- Bonfante P (2007). Laser microdissection reveals that transcripts for five plant and one fungal phosphate transporter genes are contemporaneously present in arbusculated cells. *Mol Plant Microbe Interact* 20, 1055-1062.
- Bonin S et al. (2005). DNA and RNA obtained from Bouin's fixed tissues. *J Clin Pathol* 58, 313-316.
- Bonini F et al. (2007). Surface imprinted beads for the recognition of human serum albumin. *Biosensors and Bioelectronics* 22, 2322-2328.
- Boopathi E et al. (2004). Regulation of murine cytochrome c oxidase Vb gene expression during myogenesis: yy-1 and heterogeneous nuclear ribonucleoprotein d-like protein (JKTBP1) reciprocally regulate transcription activity by physical interaction with the BERF-1/zbp-89 factor. *J Biol Chem* 279, 35242-35254.
- Bosco MC et al. (2004). Hypoxia inhibits the expression of the CCR5 chemokine receptor in macrophages. *Cell Immunol* 228, 1-7.
- Bosco MC et al. (2004). Hypoxia Selectively Inhibits Monocyte Chemoattractant Protein-1 Production by Macrophages. *J Immunol* 172, 1681-1690.
- Boulanger P (2007). The 3-O-(3-PRIME;-3-PRIME;-dimethylsuccinyl) derivative of betulinic acid (DSB) inhibits the assembly of virus-like particles in HIV-1 Gagprecursor-expressing cells. *Antiviral Ther* 12, 1185-1203.
- Bousfield GR et al. (2007). All-or-none N-glycosylation in primate follicle-stimulating hormone [beta]-subunits. *Mol Cell Endocrinol* 260-262, 40-48.
- Bowers AJ and Boylan JF (2004). Nek8, a NIMA family kinase member, is overexpressed in primary human breast tumors. *Gene* 328, 135-142.
- Boye-Harnasch M and Cullin C (2006). A novel in vitro filter trap assay identifies tannic acid as an amyloid aggregation inducer for HET-s. *J Biotech* 125, 222-230.
- Bozzo F et al. (2007). Antiproliferative effect of conjugated linoleic acid in caco-2 cells: Involvement of PPARγ and APC/β-catenin pathways. *Chemico-Biol Interact* 169, 110-121.
- Brown JM et al. (2005). Effects of roflumilast on silica-exacerbated systemic autoimmune disease in New Zealand mixed mice. *Am J Physiol Lung Cell Mol Physiol* 289, L990-L998.
- Brun P et al. (2005). Exposure to bacterial cell wall products triggers an inflammatory phenotype in hepatic stellate cells. *Am J Physiol Gastrointest Liver Physiol* 289, G571-G578.
- Brun P et al. (2007). Increased intestinal permeability in obese mice: new evidence in the pathogenesis of nonalcoholic steatohepatitis. *Am J Physiol Gastrointest Liver Physiol* 292, G518-G525.
- Brunelleschi S (2007). Minor polar compounds extra-virgin olive oil extract (MPC-OOE) inhibits NF-κB translocation in human monocyte/macrophages. *Pharmacol Res* 56, 542-549.
- Buijs JT et al. (2007). BMP7, a Putative Regulator of Epithelial Homeostasis in the Human Prostate, Is a Potent Inhibitor of Prostate Cancer Bone Metastasis in Vivo. *Am J Pathol* 171, 1047-1057.
- Bukshpan S (2007). SDS-PAGE focusing: Preparative aspects. *Anal Chem* 79, 8624-8630.
- Bunn RC et al. (2004). IGFBP-4 degradation by pregnancy-associated plasma protein-A in MC3T3 osteoblasts. *Biochem Biophys Res Commun* 325, 698-706.
- Buschmann A et al. (2004). Neuronal accumulation of abnormal prion protein in sheep carrying a scrapie-resistant genotype (PrPARR/ARR). *J Gen Virol* 85, 2727-2733.
- Buschmann A et al. (2006). Atypical BSE in Germany—Proof of transmissibility and biochemical characterization. *Vet Microbiol* 117, 103-116.
- Butler J et al. (2007). Phospholamban and sarcolipin are maintained in the endoplasmic reticulum by retrieval from the ER-Golgi intermediate compartment. *Cardiovasc Res* 74, 114-123.
- Camprostrini N et al. (2004). Proteomic analysis of an orthotopic neuroblastoma xenograft animal model. *J Chromatogr B* 808, 279-286.
- Cao A et al. (2007). Seasonal variations in haemocyte response in the mussel *Mytilus galloprovincialis* Lmk. *Aquaculture* 263, 310-319.
- Carattino MD et al. (2005). Epithelial Sodium Channel Inhibition by AMP-activated Protein Kinase in Oocytes and Polarized Renal Epithelial Cells. *J Biol Chem* 280, 17608-17616.
- Carboni L et al. (2006). Proteomic analysis of rat hippocampus after repeated psychosocial stress. *Neuroscience* 137, 1237-1246.
- Carboni L et al. (2006). Proteomic analysis of rat hippocampus and frontal cortex after chronic treatment with fluoxetine or putative novel antidepressants: CRF1 and NK1 receptor antagonists. *Eur Neuropsychopharmacol* 16, 521-537.
- Cascio P (2007). Enhanced expression of interferon-γ-induced antigen-processing machinery components in a spontaneously occurring cancer. *Neoplasia* 9, 960-969.
- Castagliuolo I et al. (2006). Mucosal genetic immunization against four adhesins protects against *Staphylococcus aureus*-induced mastitis in mice. *Vaccine* 24, 4393-4402.
- Cayuso J et al. (2006). The Sonic hedgehog pathway independently controls the patterning, proliferation and survival of neuroepithelial cells by regulating Gli activity. *Development* 133, 517-528.
- Cecconi D et al. (2007). Proteomic analysis of rat cortical neurons after fluoxetine treatment. *Brain Res* 1135, 41-51.
- Ceolotto G et al. (2006). Insulin generates free radicals in human fibroblasts ex vivo by a protein kinase C-dependent mechanism, which is inhibited by pravastatin. *Free Rad Biol Med* 41, 473-483.
- Cervený KL and Jensen RE (2003). The WD-repeats of Net2p Interact with Dnm1p and Fis1p to Regulate Division of Mitochondria. *Mol Biol Cell* 14, 4126-4139.
- Chakraborty S et al. (2006). α-Actinin 4 Potentiates Myocyte Enhancer Factor-2 Transcription Activity by Antagonizing Histone Deacetylase 7. *J Biol Chem* 281, 35070-35080.
- Cha-Molstad H et al. (2007). The interaction of C-Rel with C/EBPβ enhances C/EBPβ binding to the C-reactive protein gene promoter. *Mol Immunol* 44, 2933-2942.
- Chandra J et al. (2007). Interaction of *Candida albicans* with Adherent Human Peripheral Blood Mononuclear Cells Increases *C. albicans* Biofilm Formation and Results in Differential Expression of Pro- and Anti-Inflammatory Cytokines. *Infect Immun* 75, 2612-2620.
- Chang IY et al. (2005). Small interfering RNA-induced suppression of ERCC1 enhances sensitivity of human cancer cells to cisplatin. *Biochem Biophys Res Commun* 327, 225-233.
- Chau CH et al. (2005). Etk/Bmx mediates expression of stress-induced adaptive genes VEGF, PAI-1, and iNOS via multiple signaling cascades in different cell systems. *Am J Physiol Cell Physiol* 289, C444-C454.
- Chaudhary A et al. (2007). Differential protein expression of peroxiredoxin I and II by benzo(a)pyrene and quercetin treatment in 22Rv1 and PrEC prostate cell lines. *Toxicol Appl Pharmacol* 220, 197-210.
- Chen G et al. (2005). Modulation of Androgen Receptor Transactivation by FoxH1: A Newly Identified Androgen Receptor Corepressor. *J Biol Chem* 280, 36355-36363.

- Chen JR et al. (2005). Transient Versus Sustained Phosphorylation and Nuclear Accumulation of ERKs Underlie Anti-Versus Pro-apoptotic Effects of Estrogens. *J Biol Chem* 280, 4632-4638.
- Chen LH et al. (2007). Thapsigargin sensitizes human melanoma cells to TRAIL-induced apoptosis by up-regulation of TRAIL-R2 through the unfolded protein response. *Carcinogenesis* 28, 2328-2336.
- Chen WS and Bear MF (2007). Activity-dependent regulation of NR2B translation contributes to metaplasticity in mouse visual cortex. *Neuropharmacology* 52, 200-214.
- Cheng LT et al. (2005). Atypical Fusion Peptide of Nelson Bay Virus Fusion-Associated Small Transmembrane Protein. *J Virol* 79, 1853-1860.
- Cheroni C et al. (2005). Accumulation of human SOD1 and ubiquitinated deposits in the spinal cord of SOD1G93A mice during motor neuron disease progression correlates with a decrease of proteasome. *Neurobiol Dis* 18, 509-522.
- Chertova E et al. (2006). Proteomic and Biochemical Analysis of Purified Human Immunodeficiency Virus Type 1 Produced from Infected Monocyte-Derived Macrophages. *J Virol* 80, 9039-9052.
- Cibelli G (2007). Radiofrequency radiation (900 MHz) induces Egr-1 gene expression unaffected cell-cycle control in human neuroblastoma cells. *J Cell Physiol* 213, 759-767.
- Ciucci A et al. (2006). Induction of Apoptosis in Estrogen Receptor-Negative Breast Cancer Cells by Natural and Synthetic Cyclopentenones: Role of the I $\kappa$ B Kinase/Nuclear Factor- $\kappa$ B Pathway. *Mol Pharmacol* 70, 1812-1821.
- Clavijo C et al. (2007). Protein kinase C $\delta$ -dependent and -independent signaling in genotoxic response to treatment of desferrioxamine, a hypoxia-mimetic agent. *Am J Physiol Cell Physiol* 292, C2150-C2160.
- Claypool SM et al. (2006). Mitochondrial mislocalization and altered assembly of a cluster of Barth syndrome mutant tafazzins. *J Cell Biol* 174, 379-390.
- Clifton JG et al. (2006). Identification of members of the annexin family in the detergent-insoluble fraction of rat Morris hepatoma plasma membranes. *J Chromatogr A* 1123, 205-211.
- Coelho JE et al. (2006). Hypoxia-induced desensitization and internalization of adenosine A1 receptors in the rat hippocampus. *Neuroscience* 138, 1195-1203.
- Collier JJ et al. (2006). Pro- and Antiapoptotic Proteins Regulate Apoptosis but Do Not Protect Against Cytokine-Mediated Cytotoxicity in Rat Islets and  $\beta$ -Cell Lines. *Diabetes* 55, 1398-1406.
- Cooper J et al. (2006). Evaluation of the anterior cruciate ligament, medial collateral ligament, achilles tendon and patellar tendon as cell sources for tissue-engineered ligament. *Biomaterials* 27, 2747-2754.
- Cordier C et al. (2006). Transmission and characterization of bovine spongiform encephalopathy sources in two ovine transgenic mouse lines (TgOvPrP4 and TgOvPrP59). *J Gen Virol* 87, 3763-3771.
- Cordle A and Landreth G (2005). 3-Hydroxy-3-Methylglutaryl-Coenzyme A Reductase Inhibitors Attenuate  $\beta$ -Amyloid-Induced Microglial Inflammatory Responses. *J Neurosci* 25, 299-307.
- Cordle A et al. (2005). Mechanisms of Statin-mediated Inhibition of Small G-protein Function. *J Biol Chem* 280, 34202-34209.
- Corey DA and Kelley TJ (2007). Elevated small GTPase activation influences the cell proliferation signaling control in Niemann-Pick type C fibroblasts. *Biochimica Biophys Acta (BBA) - Mol Basis Dis* 1772, 748-754.
- Cotney J et al. (2007). Relative abundance of the human mitochondrial transcription system and distinct roles for h-mtTFB1 and h-mtTFB2 in mitochondrial biogenesis and gene expression. *Nucleic Acids Res* 35, 4042-4054.
- Crigler L et al. (2006). Human mesenchymal stem cell subpopulations express a variety of neuro-regulatory molecules and promote neuronal cell survival and neurogenesis. *Exp Neurol* 198, 54-64.
- Crusius K et al. (2006). Tab2, a novel recombinant polypeptide tag offering sensitive and specific protein detection and reliable affinity purification. *Gene* 380, 111-119.
- Cunha GMA et al. (2006). Increased density and synapto-protective effect of adenosine A2A receptors upon sub-chronic restraint stress. *Neuroscience* 141, 1775-1781.
- Cunha-Oliveira T et al. (2006). Mitochondrial dysfunction and caspase activation in rat cortical neurons treated with cocaine or amphetamine. *Brain Res* 1089, 44-54.
- Cushman I et al. (2006). Using peptide arrays to define nuclear carrier binding sites on nucleoporins. *Methods* 39, 329-341.
- Dagnon K et al. (2005). Expression of Erythropoietin and Erythropoietin Receptor in Non-Small Cell Lung Carcinomas. *Clin Cancer Res* 11, 993-999.
- Dai Q et al. (2005). Ethanol alters cellular activation and CD14 partitioning in lipid rafts. *Biochem Biophys Res Commun* 332, 37-42.
- Danzer KM et al. (2007). Different Species of  $\alpha$ -Synuclein Oligomers Induce Calcium Influx and Seeding. *J Neurosci* 27, 9220-9232.
- Darnell JC et al. (2005). Kissing complex RNAs mediate interaction between the Fragile-X mental retardation protein KH2 domain and brain polyribosomes. *Genes and Dev* 19, 903-918.
- Das S et al. (2006). Activation of the RNA-dependent Protein Kinase PKR Promoter in the Absence of Interferon Is Dependent Upon Sp Proteins. *J Biol Chem* 281, 3244-3253.
- Das SC and Pattnaik AK (2004). Phosphorylation of Vesicular Stomatitis Virus Phosphoprotein P Is Indispensable for Virus Growth. *J Virol* 78, 6420-6430.
- Das SC and Pattnaik AK (2005). Role of the Hypervariable Hinge Region of Phosphoprotein P of Vesicular Stomatitis Virus in Viral RNA Synthesis and Assembly of Infectious Virus Particles. *J Virol* 79, 8101-8112.
- Davies RA et al. (2006). Hepatic oval cell response to the choline-deficient, ethionine supplemented model of murine liver injury is attenuated by the administration of a cyclo-oxygenase 2 inhibitor. *Carcinogenesis* 27, 1607-1616.
- Davis J et al. (2004). Early-onset and Robust Cerebral Microvascular Accumulation of Amyloid  $\beta$ -Protein in Transgenic Mice Expressing Low Levels of a Vasculotropic Dutch/Iowa Mutant Form of Amyloid  $\beta$ -Protein Precursor. *J Biol Chem* 279, 20296-20306.
- Davis J et al. (2006). Deficient cerebral clearance of vasculotropic mutant Dutch/Iowa Double A[ss] in human A[ss]PP transgenic mice. *Neurobiol Aging* 27, 946-954.
- Davis MR et al. (2004). Thrombus-induced endothelial dysfunction: Hemoglobin and fibrin decrease nitric oxide bioactivity without altering eNOS. *J Surg Res* 122, 121-129.
- Demoz M et al. (2006). High yield synthesis and characterization of phosphorylated recombinant human procathepsin D expressed in mammalian cells. *Protein Expression and Purification* 45, 157-167.
- Denkins Y et al. (2005). Role of  $\omega$ -3 polyunsaturated fatty acids on cyclooxygenase-2 metabolism in brain-metastatic melanoma. *J Lipid Res* 46, 1278-1284.
- Dey A et al. (2005). Evidence for cytochrome P450 2E1 catalytic activity and expression in rat blood lymphocytes. *Life Sci* 77, 1082-1093.
- Dey A et al. (2006). Evidence for cytochrome P450 3A expression and catalytic activity in rat blood lymphocytes. *Life Sci* 79, 175-183.
- Dong X et al. (2004). In situ dynamically monitoring the proteolytic function of the ubiquitin-proteasome system in cultured cardiac myocytes. *Am J Physiol Heart Circ Physiol* 287, H1417-H1425.
- Doyle J et al. (2006). Two Domains That Control Prefusion Stability and Transport Competence of the Measles Virus Fusion Protein. *J Virol* 80, 1524-1536.
- Du L et al. (2004). Inhibition of Cell Proliferation and Cell Cycle Progression by Specific Inhibition of Basal JNK Activity: Evidence That Mitotic Bcl-2 Phosphorylation Is Jnk-Independent. *J Biol Chem* 279, 11957-11966.
- Duarte AI et al. (2005). Insulin neuroprotection against oxidative stress in cortical neurons--Involvement of uric acid and glutathione antioxidant defenses. *Free Rad Biol Med* 39, 876-889.
- Duarte J (2007). Modification of purinergic signaling in the hippocampus of streptozotocin-induced diabetic rats. *Neuroscience* 149, 382-391.
- Duarte JMN et al. (2006). Modification of adenosine A1 and A2A receptor density in the hippocampus of streptozotocin-induced diabetic rats. *Neurochem Int* 48, 144-150.
- Duarte JMN et al. (2007). Increase of cannabinoid CB1 receptor density in the hippocampus of streptozotocin-induced diabetic rats. *Exp Neurol* 204, 479-484.

- Dunn CD et al. (2006). A Genomewide Screen for Petite-negative Yeast Strains Yields a New Subunit of the i-AAA Protease Complex. *Mol Biol Cell* 17, 213-226.
- Durand SH et al. (2006). Lipoteichoic Acid Increases TLR and Functional Chemokine Expression while Reducing Dentin Formation in In Vitro Differentiated Human Odontoblasts. *J Immunol* 176, 2880-2887.
- Efanov AM et al. (2005). A Novel Glucokinase Activator Modulates Pancreatic Islet and Hepatocyte Function. *Endocrinology* 146, 3696-3701.
- Eiden M et al. (2006). Synergistic and strain-specific effects of bovine spongiform encephalopathy and scrapie prions in the cell-free conversion of recombinant prion protein. *J Gen Virol* 87, 3753-3761.
- El Karkouri K et al. (2007). Identification of Internal Transcribed Spacer Sequence Motifs in Truffles: a First Step toward Their DNA Bar Coding. *Appl Environ Microbiol* 73, 5320-5330.
- Fahmy K et al. (2006). Secondary Structure and Pd(II) Coordination in S-Layer Proteins from *Bacillus sphaericus* Studied by Infrared and X-Ray Absorption Spectroscopy. *Biophys J* 91, 996-1007.
- Fairn GD and McMaster CR (2005). Studying phospholipid metabolism using yeast systematic and chemical genetics. *Methods* 36, 102-108.
- Fairn GD et al. (2007). A Chemogenomic Screen in *Saccharomyces cerevisiae* Uncovers a Primary Role for the Mitochondria in Farnesol Toxicity and Its Regulation by the Pkc1 Pathway. *J Biol Chem* 282, 4868-4874.
- Fan Q et al. (2004). Candesartan reduced advanced glycation end-products accumulation and diminished nitro-oxidative stress in type 2 diabetic KK/Ta mice. *Nephrol Dial Transplant* 19, 3012-3020.
- Fan R et al. (2007). Minocycline Reduces Microglial Activation and Improves Behavioral Deficits in a Transgenic Model of Cerebral Microvascular Amyloid. *J Neurosci* 27, 3057-3063.
- Fang J et al. (2007). Characterization of Chimeric *Bacillus thuringiensis* Vip3 Toxins. *Appl Environ Microbiol* 73, 956-961.
- Farley SM et al. (2006). Menin Localizes to Chromatin Through an ATR-CHK1 Mediated Pathway After UV-Induced DNA Damage. *J Surg Res* 133, 29-37.
- Ferrari F et al. (2007). The fragile X mental retardation protein-RNP granules show an mGluR-dependent localization in the post-synaptic spines. *Mol Cell Neurosci* 34, 343-354.
- Ferreiro E et al. (2006). An endoplasmic-reticulum-specific apoptotic pathway is involved in prion and amyloid-beta peptides neurotoxicity. *Neurobiol Dis* 23, 669-678.
- Fettucciari K et al. (2006). Group B *Streptococcus* Induces Macrophage Apoptosis by Calpain Activation. *J Immunol* 176, 7542-7556.
- Fifre A et al. (2006). Microtubule-associated Protein MAP1A, MAP1B, and MAP2 Proteolysis during Soluble Amyloid  $\beta$ -Peptide-induced Neuronal Apoptosis: Synergistic Involvement Of Calpain And Caspase-3. *J Biol Chem* 281, 229-240.
- Filigheddu N et al. (2007). Ghrelin and Des-Acyl Ghrelin Promote Differentiation and Fusion of C2C12 Skeletal Muscle Cells. *Mol Biol Cell* 18, 986-994.
- Filipkowski P et al. (2007). Identification, cloning, expression, and characterization of a highly thermostable single-stranded-DNA-binding protein (SSB) from *Deinococcus murrayi*. *Protein Expression and Purification* 53, 201-208.
- Fleury C et al. (2006). Effect of cobalt and chromium ions on human MG-63 osteoblasts in vitro: Morphology, cytotoxicity, and oxidative stress. *Biomaterials* 27, 3351-3360.
- Fogdell-Hahn A (2007). Purification of infectious human herpesvirus 6A virions and association of host cell proteins. *Virology Journal* 4, U1-U11.
- Fowlkes JL et al. (2004). Regulation of Insulin-Like Growth Factor (IGF)-I Action by Matrix Metalloproteinase-3 Involves Selective Disruption of IGF-I/IGF-Binding Protein-3 Complexes. *Endocrinology* 145, 620-626.
- Fueger PT et al. (2007). Glucose kinetics and exercise tolerance in mice lacking the GLUT4 glucose transporter. *J Physiol* 582, 801-812.
- Fueger PT et al. (2007). Phosphorylation Barriers to Skeletal and Cardiac Muscle Glucose Uptakes in High-Fat Fed Mice: Studies in Mice With a 50% Reduction of Hexokinase II. *Diabetes* 56, 2476-2484.
- Fukumoto S et al. (2003). High-Level Expression of Truncated Surface Antigen P50 of *Babesia gibsoni* in Insect Cells by Baculovirus and Evaluation of Its Immunogenicity and Antigenicity. *Clin Diagn Lab Immunol* 10, 596-601.
- Galeotti N et al. (2006). Signaling pathway of morphine induced acute thermal hyperalgesia in mice. *Pain* 123, 294-305.
- Gambacorta A (2007). *Halomonas sinaiensis* sp. nov., a novel halophilic bacterium isolated from a salt lake inside Ras Muhammad Park, Egypt. *Extremophiles* 11, 789-796.
- Gambarana C (2007). Behavioral expression of cocaine sensitization in rats is accompanied by a distinct pattern of modifications in the PKA/DARPP-32 signaling pathway. *J Neurochem* 103, 1168-1183.
- Gao X et al. (2007). TNF- $\alpha$  Contributes to Endothelial Dysfunction by Upregulating Arginase in Ischemia/Reperfusion Injury. *Arterioscler Thromb Vasc Biol* 27, 1269-1275.
- Gao X et al. (2007). Tumor Necrosis Factor- $\alpha$  Induces Endothelial Dysfunction in Leprdb Mice. *Circulation* 115, 245-254.
- Garcia RC et al. (2005). Complement Receptor 3 Binds the *Borrelia burgdorferi* Outer Surface Proteins OspA and OspB in an iC3b-Independent Manner. *Infect Immun* 73, 6138-6142.
- George CX et al. (2005). Expression of Interferon-inducible RNA Adenosine Deaminase ADAR1 during Pathogen Infection and Mouse Embryo Development Involves Tissue-selective Promoter Utilization and Alternative Splicing. *J Biol Chem* 280, 15020-15028.
- Gharavi NM et al. (2006). Role of Endothelial Nitric Oxide Synthase in the Regulation of SREBP Activation by Oxidized Phospholipids. *Circ Res* 98, 768-776.
- Gil J et al. (2003). Cytosolic and mitochondrial ROS in staurosporine-induced retinal cell apoptosis. *Free Rad Biol Med* 35, 1500-1514.
- Gillespie AS and Elliott E (2005). Comparative advantages of imidazole-sodium dodecyl sulfate-zinc reverse staining in polyacrylamide gels. *Analytical Biochem* 345, 158-160.
- Giunti R et al. (2007). Both translocon and a cation channel are involved in the passive Ca<sup>2+</sup> leak from the endoplasmic reticulum: A mechanistic study on rat liver microsomes. *Arch Biochem Biophys* 462, 115-121.
- Giuriso E et al. (2003). T Cell Receptor Can Be Recruited to a Subset of Plasma Membrane Rafts, Independently of Cell Signaling and Attending to Raft Clustering. *J Biol Chem* 278, 6771-6778.
- Glatzel M et al. (2003). Extraneural Pathologic Prion Protein in Sporadic Creutzfeldt-Jakob Disease. *N Engl J Med* 349, 1812-1820.
- Goda S et al. (2006). Matrix Metalloproteinase-1 Produced by Human CXCL12-Stimulated Natural Killer Cells. *Am J Pathol* 169, 445-458.
- Gomelsky L et al. (2003). Identification and in vivo characterization of PpaA, a regulator of photosystem formation in *Rhodobacter sphaeroides*. *Microbiology* 149, 377-388.
- Gonzalez F (2007). P2Y(2) nucleotide receptors inhibit trauma-induced death of astrocytic cells. *J Neurochem* 103, 1785-1800.
- Gonzalez-Riopadre M et al. (2007). Effect of thermal stress on protein expression in the mussel *Mytilus galloprovincialis* Lmk. *Comp Biochem Physiol Part B: Biochem Mol Biol* 147, 531-540.
- Govers R et al. (2004). Insulin Increases Cell Surface GLUT4 Levels by Dose Dependently Discharging GLUT4 into a Cell Surface Recycling Pathway. *Mol Cell Biol* 24, 6456-6466.
- Graat HCA et al. (2007). Enhanced tumor cell kill by combined treatment with a small-molecule antagonist of mouse double minute 2 and adenoviruses encoding p53. *Mol Cancer Ther* 6, 1552-1561.
- Graham NA and Asthagiri AR (2004). Epidermal Growth Factor-mediated T-cell Factor/Lymphoid Enhancer Factor Transcriptional Activity Is Essential but Not Sufficient for Cell Cycle Progression in Nontransformed Mammary Epithelial Cells. *J Biol Chem* 279, 23517-23524.
- Graham RI et al. (2004). Genetically variable nucleopolyhedroviruses isolated from spatially separate populations of the winter moth *Operophtera brumata* (Lepidoptera: Geometridae) in Orkney. *J Invertebr Pathol* 87, 29-38.
- Grande F et al. (2007). Synthesis and antitumor activities of a series of novel quinoxalinhydrazides. *Bioorganic & Med Chem* 15, 288-294.



- Graybill ER et al. (2007). Functional comparison of citrate synthase isoforms from *S. cerevisiae*. *Arch Biochem Biophys* 465, 26-37.
- Grontvedt RN and Espelid S (2004). Vaccination and immune responses against atypical *Aeromonas salmonicida* in spotted wolffish (*Anarhichas minor Olafsen*) juveniles. *Fish & Shellfish Immunol* 16, 271-285.
- Gu ML et al. (2003). Development of an animal-human antibody complex for use as a control in ELISA. *J Pharm Biomed Analysis* 32, 523-529.
- Guarna M et al. (2003). Anti-mu opioid antiserum against the third external loop of the cloned mu-opioid receptor acts as a mu receptor neutral antagonist. *Mol Brain Res* 119, 100-110.
- Guilbault B and Kay RJ (2004). RasGRP1 Sensitizes an Immature B Cell Line to Antigen Receptor-induced Apoptosis. *J Biol Chem* 279, 19523-19530.
- Gupta SC et al. (2007). Adverse effect of organophosphate compounds, dichlorvos and chlorpyrifos in the reproductive tissues of transgenic *Drosophila melanogaster*: 70 kDa heat shock protein as a marker of cellular damage. *Toxicology* 238, 1-14.
- Gurevich V (2007). Cone arrestin binding to JNK3 and Mdm2: conformational preference and localization of interaction sites. *J Neurochem* 103, 1053-1062.
- Guzzi A (2007). Stromal complement receptor CD21/35 facilitates lymphoid prion colonization and pathogenesis. *J Immunol* 179, 6144-6152.
- Hachiya NS et al. (2005). More than a 100-fold increase in immunoblot signals of laser-microdissected inclusion bodies with an excessive aggregation property by oligomeric actin interacting protein 2/d-lactate dehydrogenase protein 2. *Analytical Biochem* 347, 106-111.
- Haitcock E et al. (2005). From the Cover: Age-related changes of nuclear architecture in *Caenorhabditis elegans*. *Proc Natl Acad Sci USA* 102, 16690-16695.
- Hall A (2007). Activin acts with nerve growth factor to regulate calcitonin gene-related peptide mRNA in sensory neurons. *Neuroscience* 150, 665-674.
- Hallows KR et al. (2006). Up-regulation of AMP-activated Kinase by Dysfunctional Cystic Fibrosis Transmembrane Conductance Regulator in Cystic Fibrosis Airway Epithelial Cells Mitigates Excessive Inflammation. *J Biol Chem* 281, 4231-4241.
- Handy DE et al. (2006). Aminoglycosides Decrease Glutathione Peroxidase-1 Activity by Interfering with Selenocysteine Incorporation. *J Biol Chem* 281, 3382-3388.
- Hanson SM et al. (2007). Each rhodopsin molecule binds its own arrestin. *Proc Natl Acad Sci USA* 104, 3125-3128.
- Harrison-Bernard LM et al. (2006). Efferent arterioles exclusively express the subtype 1A angiotensin receptor: functional insights from genetic mouse models. *Am J Physiol Renal Physiol* 290, F1177-F1186.
- Henning P et al. (2006). Adenovirus type 5 fiber knob domain has a critical role in fiber protein synthesis and encapsidation. *J Gen Virol* 87, 3151-3160.
- Hersey P (2007). Inhibition of MEK sensitizes human melanoma cells to endoplasmic reticulum stress-induced apoptosis. *Cancer Res* 67, 9750-9761.
- Hersey P (2007). Regulation of docetaxel-induced apoptosis of human melanoma cells by different isoforms of protein kinase C. *Mol Cancer Res* 5, 1073-1081.
- Hersey P (2007). Temozolomide induces senescence but not apoptosis in human melanoma cells. *Br J Cancer* 97, 1225-1233.
- Hicks SW and Machamer CE (2005). Isoform-specific interaction of Golgin-160 with the Golgi-associated protein PIST. *J Biol Chem* 280, 28944-28951.
- Higashibata A et al. (2006). Decreased expression of myogenic transcription factors and myosin heavy chains in *Caenorhabditis elegans* muscles developed during spaceflight. *J Exp Biol* 209, 3209-3218.
- Hirumura M et al. (2006). Identification of Nerve Growth Factor-responsive Element of the TCL1 Promoter as a Novel Negative Regulatory Element. *J Biol Chem* 281, 27753-27764.
- Ho YP et al. (2006). Evaluating the intracellular stability and unpacking of DNA nanocomplexes by quantum dots-FRET. *J Controlled Release* 116, 83-89.
- Hofmann G (2007). Is cold the new hot? Elevated ubiquitin-conjugated protein levels in tissues of Antarctic fish as evidence for cold-denaturation of proteins in vivo. *J Comp Physiol B-Biochem Systemic Environ Physiol* 177, 857-866.
- Hong SS et al. (2003). Identification of Adenovirus (Ad) Penton Base Neutralizing Epitopes by Use of Sera from Patients Who Had Received Conditionally Replicative Ad (Add1520) for Treatment of Liver Tumors. *J Virol* 77, 10366-10375.
- Hoos MD et al. (2007). Inhibition of Familial Cerebral Amyloid Angiopathy Mutant Amyloid beta-Protein Fibril Assembly by Myelin Basic Protein. *J Biol Chem* 282, 9952-9961.
- Hsu YH et al. (2006). Sumoylated SnoN Represses Transcription in a Promoter-specific Manner. *J Biol Chem* 281, 33008-33018.
- Hua NP et al. (2007). *Halobacillus profundus* sp. nov. and *Halobacillus kuroshimensis* sp. nov., moderately halophilic bacteria isolated from a deep-sea methane cold seep. *Int J Syst Evol Microbiol* 57, 1243-1249.
- Huang W et al. (2003). A broad-spectrum peptide inhibitor of  $\beta$ -lactamase identified using phage display and peptide arrays. *Protein Eng* 16, 853-860.
- Hurtado CAR and Rachubinski RA (2002). Isolation and Characterization of YIBEM1, a Gene Required for Cell Polarization and Differentiation in the Dimorphic Yeast *Yarrowia lipolytica*. *Eukaryot Cell* 1, 526-537.
- Hurtado CAR and Rachubinski RA (2002). YIBMH1 encodes a 14-3-3 protein that promotes filamentous growth in the dimorphic yeast *Yarrowia lipolytica*. *Microbiology* 148, 3725-3735.
- Hwang I (2007). Effect of fiber type on postmortem proteolysis in longissimus muscle of Landrace and Korean native black pigs. *Meat Sci* 77, 482-491.
- Israfi DA et al. (2007). Cardamonin inhibits COX and iNOS expression via inhibition of p65NF- $\kappa$ B nuclear translocation and  $\kappa$ B phosphorylation in RAW 264.7 macrophage cells. *Mol Immunol* 44, 673-679.
- Iltner LM et al. (2004). Aspartate69 of the calcitonin-like receptor is required for its functional expression together with receptor-activity-modifying proteins 1 and -2. *Biochem Biophys Res Commun* 319, 1203-1209.
- James ER et al. (2003). Presence of a Transcriptionally Active Glucocorticoid Receptor  $\alpha$  in Lens Epithelial Cells. *Invest Ophthalmol Vis Sci* 44, 5269-5276.
- Jang-Lee J et al. (2007). Glycomics Analysis of *Schistosoma mansoni* Egg and Cercarial Secretions. *Mol Cell Proteomics* 6, 1485-1499.
- Jenny MJ et al. (2006). Regulation of metallothionein genes in the American oyster (*Crassostrea virginica*): Ontogeny and differential expression in response to different stressors. *Gene* 379, 156-165.
- Jensen MV et al. (2006). Compensatory Responses to Pyruvate Carboxylase Suppression in Islet beta-Cells: Preservation Of Glucose-Stimulated Insulin Secretion. *J Biol Chem* 281, 22342-22351.
- Jeyaraju DV et al. (2006). Phosphorylation and cleavage of presenilin-associated rhomboid-like protein (PARL) promotes changes in mitochondrial morphology. *Proc Natl Acad Sci USA* 103, 18562-18567.
- Jia D et al. (2006). Glucocorticoids Act Directly on Osteoclasts to Increase Their Life Span and Reduce Bone Density. *Endocrinology* 147, 5592-5599.
- Jiang CC et al. (2007). Tunicamycin Sensitizes Human Melanoma Cells to Tumor Necrosis Factor-Related Apoptosis-Inducing Ligand-Induced Apoptosis by Up-regulation of TRAIL-R2 via the Unfolded Protein Response. *Cancer Res* 67, 5880-5888.
- Johnson BM et al. (2006). Characterization Of Transport Protein Expression In Multidrug Resistance-Associated Protein (MRP) 2-Deficient Rats. *Drug Metab Dispos* 34, 556-562.
- Johri A et al. (2006). Effect of prenatal exposure of deltamethrin on the ontogeny of xenobiotic metabolizing cytochrome P450s in the brain and liver of offspring. *Toxicol Appl Pharmacol* 214, 279-289.
- Johri A et al. (2006). Long lasting effects of prenatal exposure to deltamethrin on cerebral and hepatic cytochrome P450s and behavioral activity in rat offspring. *Eur J Pharmacol* 544, 58-68.
- Kalra N et al. (2007). Theaflavins induced apoptosis of LNCaP cells is mediated through induction of p53, down-regulation of NF- $\kappa$ B and mitogen-activated protein kinases pathways. *Life Sci* 80, 2137-2146.
- Kang X et al. (2007). Adipogenesis of murine embryonic stem cells in a three-dimensional culture system using electrospun polymer scaffolds. *Biomaterials* 28, 450-458.
- Kapoor N et al. (2006). Cytochrome P450 1A isoenzymes in brain cells: Expression and inducibility in cultured rat brain neuronal and glial cells. *Life Sci* 79, 2387-2394.

- Kapoor N et al. (2006). Differences in sensitivity of cultured rat brain neuronal and glial cytochrome P450 2E1 to ethanol. *Life Sci* 79, 1514-1522.
- Kashlan OB et al. (2007). Small Heat Shock Protein  $\alpha$ A-crystallin Regulates Epithelial Sodium Channel Expression. *J Biol Chem* 282, 28149-28156.
- Kato S (2007). Increased susceptibility of small intestine to NSAID-provoked ulceration in rats with adjuvant-induced arthritis: Involvement of enhanced expression of TLR4. *Life Sci* 81, 1309-1316.
- Kawai S et al. (2006). Odd-Skipped Related 2 gene transcription is regulated by CCAAT enhancer-binding protein  $\delta$  in mesenchymal C3H10T1/2 cells. *Genes Cells* 11, 163-175.
- Kawasaki S et al. (2006). Clusters of Corneal Epithelial Cells Reside Ectopically in Human Conjunctival Epithelium. *Invest Ophthalmol Vis Sci* 47, 1359-1367.
- Kawasaki S et al. (2006). Expression and tissue distribution of p63 isoforms in human ocular surface epithelia. *Exp Eye Res* 82, 293-299.
- Kawate H et al. (2005). Impaired Nuclear Translocation, Nuclear Matrix Targeting, and Intracellular Mobility of Mutant Androgen Receptors Carrying Amino Acid Substitutions in the Deoxyribonucleic Acid-Binding Domain Derived from Androgen Insensitivity Syndrome Patients. *J Clin Endocrinol Metab* 90, 6162-6169.
- Kawate H et al. (2007). Mutual transactivational repression of Runx2 and the androgen receptor by an impairment of their normal compartmentalization. *J Steroid Biochem Mol Biol* 105, 46-56.
- Kepler CJ et al. (2005). Sublethal effects of the toxic alga *Heterosigma akashiwo* on the southeastern oyster (*Crassostrea virginica*). *Harmful Algae* 4, 275-285.
- Kern ER et al. (2005). In Vitro Activity and Mechanism of Action of Methylenecyclopropane Analogs of Nucleosides against Herpesvirus Replication. *Antimicrob Agents Chemother* 49, 1039-1045.
- Kessler-Icekson G (2007). The Popdc gene family in the rat: Molecular cloning, characterization and expression analysis in the heart and cultured cardiomyocytes. *BIOCHIMICA BIOPHYS ACTA-GENE STRUCTURE AND EXPRESSION* 1769, 586-592.
- Kessler-Icekson G et al. (2006). Expression of procollagen C-proteinase enhancer-1 in the remodeling rat heart is stimulated by aldosterone. *Int J Biochem Cell Biol* 38, 358-365.
- Khuda S et al. (2004). The Sac3 Homologue SHD1 Is Involved in Mitotic Progression in Mammalian Cells. *J Biol Chem* 279, 46182-46190.
- Kim DJ et al. (2006). Homocysteine enhances apoptosis in human bone marrow stromal cells. *Bone* 39, 582-590.
- Kim E et al. (2003). Disruption of TR4 Orphan Nuclear Receptor Reduces the Expression of Liver Apolipoprotein E/C-I/C-II Gene Cluster. *J Biol Chem* 278, 46919-46926.
- Kim KY et al. (2007). Molecular characterization and mRNA expression during metal exposure and thermal stress of copper/zinc- and manganese-superoxide dismutases in disk abalone, *Haliotis discus discus*. *Fish & Shellfish Immunol* 23, 1043-1059.
- Kim M et al. (2007). Decreased Catalytic Activity of the Insulin-degrading Enzyme in Chromosome 10-Linked Alzheimer Disease Families. *J Biol Chem* 282, 7825-7832.
- Kima PE and Dunn W (2005). Exploiting calnexin expression on phagosomes to isolate *Leishmania* parasitophorous vacuoles. *Microbial Pathogenesis* 38, 139-145.
- Kimura M et al. (2007). A Critical Role for Allograft Inflammatory Factor-1 in the Pathogenesis of Rheumatoid Arthritis. *J Immunol* 178, 3316-3322.
- Kines KJ et al. (2006). Transduction of *Schistosoma mansoni* by vesicular stomatitis virus glycoprotein-pseudotyped Moloney murine leukemia retrovirus. *Exp Parasitology* 112, 209-220.
- Kinlough CL et al. (2004). MUC1 Membrane Trafficking Is Modulated by Multiple Interactions. *J Biol Chem* 279, 53071-53077.
- Kinlough CL et al. (2006). Recycling of MUC1 Is Dependent on Its Palmitoylation. *J Biol Chem* 281, 12112-12122.
- Kinney RM et al. (2006). Avian virulence and thermostable replication of the North American strain of West Nile virus. *J Gen Virol* 87, 3611-3622.
- Kinugasa Y et al. (2004). Neuroglycan C, a novel member of the neuregulin family. *Biochem Biophys Res Commun* 321, 1045-1049.
- Kjaerulf S and Jensen MR (2005). Comparison of different signal peptides for secretion of heterologous proteins in fission yeast. *Biochem Biophys Res Commun* 336, 974-982.
- Kjaerulf S et al. (2005). Alternative protein secretion: The Mam1 ABC transporter supports secretion of M-factor linked GFP in fission yeast. *Biochem Biophys Res Commun* 338, 1853-1859.
- Klungland H (2007). Altered expression of PGK1 in a family with phosphoglycerate kinase deficiency. *Muscle & Nerve* 36, 679-684.
- Koehler C (2007). A role for cytochrome c and cytochrome c peroxidase in electron shuttling from Erv1. *EMBO J* 26, 4801-4811.
- Koh JM et al. (2005).  $\alpha$ -Lipoic acid suppresses osteoclastogenesis despite increasing the receptor activator of nuclear factor  $\kappa$ B ligand/osteoprotegerin ratio in human bone marrow stromal cells. *J Endocrinol* 185, 401-413.
- Koller D et al. (2004). Selective Inactivation of Adrenomedullin over Calcitonin Gene-related Peptide Receptor Function by the Deletion of Amino Acids 14-20 of the Mouse Calcitonin-like Receptor. *J Biol Chem* 279, 20387-20391.
- Koshida R et al. (2005). Role of Focal Adhesion Kinase in Flow-Induced Dilatation of Coronary Arterioles. *Arterioscler Thromb Vasc Biol* 25, 2548-2553.
- Kosinski M et al. (2005). *C. elegans* sperm bud vesicles to deliver a meiotic maturation signal to distant oocytes. *Development* 132, 3357-3369.
- Kumar S et al. (2004). Expression of *Babesia equi* EMA-1 and EMA-2 during merozoite developmental stages in erythrocyte and their interaction with erythrocytic membrane skeleton. *Mol Biochem Parasitol* 133, 221-227.
- Kur J (2007). ADSRRS-fingerprinting and PCR MP techniques for studies of intraspecies genetic relatedness in *Staphylococcus aureus*. *J Microbiol Methods* 71, 114-122.
- Kurlawalla-Martinez C et al. (2005). Insulin Hypersensitivity and Resistance to Streptozotocin-Induced Diabetes in Mice Lacking PTEN in Adipose Tissue. *Mol Cell Biol* 25, 2498-2510.
- Kuzman JA et al. (2005). Effects of induced hyperthyroidism in normal and cardiomyopathic hamsters. *J Appl Physiol* 99, 1428-1433.
- Kuzman JA et al. (2005). L-Thyroxine activates Akt signaling in the heart. *J Mol Cell Cardiol* 39, 251-258.
- Kuzman JA et al. (2007). Rapamycin Prevents Thyroid Hormone-Induced Cardiac Hypertrophy. *Endocrinology* 148, 3477-3484.
- Kwak MK et al. (2003). Modulation of Gene Expression by Cancer Chemo-preventive Dithiolethiones through the Keap1-Nrf2 Pathway. Identification of Novel Gene Clusters For Cell Survival. *J Biol Chem* 278, 8135-8145.
- Kwak MK et al. (2007). Role of increased expression of the proteasome in the protective effects of sulforaphane against hydrogen peroxide-mediated cytotoxicity in murine neuroblastoma cells. *Free Rad Biol Med* 43, 809-817.
- Kwak MK et al. (2007). Tissue specific increase of the catalytic subunits of the 26S proteasome by indirect antioxidant dithiolethione in mice: Enhanced activity for degradation of abnormal protein. *Life Sci* 80, 2411-2420.
- Kwong WY et al. (2006). Imprinted gene expression in the rat embryo-fetal axis is altered in response to periconceptual maternal low protein diet. *Reproduction* 132, 265-277.
- Lagos D et al. (2007). Kaposi sarcoma herpesvirus-encoded vFLIP and vIRF1 regulate antigen presentation in lymphatic endothelial cells. *Blood* 109, 1550-1558.
- Lagrange M et al. (2007). Intracellular scFvs against the viral E6 oncoprotein provoke apoptosis in human papillomavirus-positive cancer cells. *Biochem Biophys Res Commun* 361, 487-492.
- Lake SL et al. (2006). Effects of body condition score at parturition and post-partum supplemental fat on metabolite and hormone concentrations of beef cows and their suckling calves. *J Anim Sci* 84, 1038-1047.
- Lamartiniere C (2007). Resveratrol suppresses prostate cancer progression in transgenic mice. *Carcinogenesis* 28, 1946-1953.
- Larance M et al. (2005). Characterization of the Role of the Rab GTPase-activating Protein AS160 in Insulin-regulated GLUT4 Trafficking. *J Biol Chem* 280, 37803-37813.

- Lareu RR et al. (2007). Collagen matrix deposition is dramatically enhanced in vitro when crowded with charged macromolecules: The biological relevance of the excluded volume effect. *FEBS Lett* 581, 2709-2714.
- Larijani M et al. (2005). Lack of MSH2 involvement differentiates V(D)J recombination from other non-homologous end joining events. *Nucleic Acids Res* 33, 6733-6742.
- Laurie NA et al. (2005). Carcinoembryonic Antigen-Related Cell Adhesion Molecule 1a-4L Suppression of Rat Hepatocellular Carcinomas. *Cancer Res* 65, 11010-11017.
- Lee JK et al. (2007). Reversible Inhibition of the Fusion Activity of Measles Virus F Protein by an Engineered Intersubunit Disulfide Bridge. *J Virol* 81, 8821-8826.
- Lee MJ et al. (2007). Feeding and Insulin Increase Leptin Translation: Importance of the Leptin Mrna Untranslated Regions. *J Biol Chem* 282, 72-80.
- Lenard NR et al. (2006). Differential coupling of beta3A- and beta3B-adrenergic receptors to endogenous and chimeric G $\alpha$ s and G $\alpha$ i. *Am J Physiol Endocrinol Metab* 291, E704-E715.
- Li AQ et al. (2007). Proteomics of the flesh fly brain reveals an abundance of upregulated heat shock proteins during pupal diapause. *J Insect Physiol* 53, 385-391.
- Li R et al. (2007). OKL38 is an oxidative stress response gene stimulated by oxidized phospholipids. *J Lipid Res* 48, 709-715.
- Li Y et al. (2006). Inducible Resistance of Tumor Cells to Tumor Necrosis Factor-Related Apoptosis-Inducing Ligand Receptor 2-Mediated Apoptosis by Generation of a Blockade at the Death Domain Function. *Cancer Res* 66, 8520-8528.
- Liang Q et al. (2006). P38 MAP Kinase Activity Is Correlated With Angiotensin II Type 1 Receptor Blocker-Induced Left Ventricular Reverse Remodeling in Spontaneously Hypertensive Heart Failure Rats. *J Cardiac Failure* 12, 479-486.
- Liedtke CM et al. (2004). Role of a PDZ1 domain of NHERF1 in the binding of airway epithelial RACK1 to NHERF1. *Am J Physiol Cell Physiol* 286, C1037-C1044.
- Lih CJ et al. (2006). Txr1: a transcriptional regulator of thrombospondin-1 that modulates cellular sensitivity to taxanes. *Genes and Dev* 20, 2082-2095.
- Lim CJ and Shen WC (2005). Comparison of monomeric and oligomeric transferrin as potential carrier in oral delivery of protein drugs. *J Controlled Release* 106, 273-286.
- Lin HY et al. (2006). Integrin  $\alpha$ v $\beta$ 3 contains a receptor site for resveratrol. *FASEB J* 20, 1742-1744.
- Liu J et al. (2005). Impairment of the ubiquitin-proteasome system in desminopathy mouse hearts. *FASEB J* 05-4869fje.
- Liu WF et al. (2006). E-cadherin engagement stimulates proliferation via Rac1. *J Cell Biol* 173, 431-441.
- Liu WF et al. (2007). Cadherins, RhoA, and Rac1 Are Differentially Required for Stretch-Mediated Proliferation in Endothelial Versus Smooth Muscle Cells. *Circ Res* 101, e44-e52.
- Louder MK et al. (2005). HIV-1 envelope pseudotyped viral vectors and infectious molecular clones expressing the same envelope glycoprotein have a similar neutralization phenotype, but culture in peripheral blood mononuclear cells is associated with decreased neutralization sensitivity. *Virology* 339, 226-238.
- Lu Q et al. (2004). EST-based genome-wide gene inactivation identifies ARAP3 as a host protein affecting cellular susceptibility to anthrax toxin. *Proc Natl Acad Sci USA* 101, 17246-17251.
- Luchetti S et al. (2006). Comparative non-radioactive RT-PCR assay: An approach to study the neurosteroids biosynthetic pathway in humans. *J Neurosci Methods* 153, 290-298.
- Lum R et al. (2004). Evidence for an Unfolding/Threading Mechanism for Protein Disaggregation by *Saccharomyces cerevisiae* Hsp104. *J Biol Chem* 279, 29139-29146.
- Lund SG et al. (2006). Turning up the heat: The effects of thermal acclimation on the kinetics of hsp70 gene expression in the eurythermal goby, *Gillichthys mirabilis*. *Comp Biochem Physiol - Part A: Mol & Integrative Physiol* 143, 435-446.
- Luo L et al. (2005). Effect of cobalt and chromium ions on MMP-1, TIMP-1, and TNF- $\alpha$  gene expression in human U937 macrophages: A role for tyrosine kinases. *Biomaterials* 26, 5587-5593.
- Maag RS et al. (2005). Caspase-resistant Golgin-160 Disrupts Apoptosis Induced by Secretory Pathway Stress and Ligation of Death Receptors. *Mol Biol Cell* 16, 3019-3027.
- Macedo T (2007). Methamphetamine induces alterations on hippocampal NMDA and AMPA receptor subunit levels and impairs spatial working memory. *Neuroscience* 150, 433-441.
- Madan M et al. (2007). Doxycycline affects diet- and bacteria-associated atherosclerosis in an ApoE heterozygote murine model: Cytokine profiling implications. *Atherosclerosis* 190, 62-72.
- Maenaka K (2007). Crystal structure of measles virus hemagglutinin provides insight into effective vaccines. *Proc Natl Acad Sci USA* 104, 19535-19540.
- Manganelli R (2007). PE is a functional domain responsible for protein translocation and localization on mycobacterial cell wall. *Mol Microbiol* 66, 1536-1547.
- Marbois B et al. (2005). Coq3 and Coq4 Define a Polypeptide Complex in Yeast Mitochondria for the Biosynthesis of Coenzyme Q. *J Biol Chem* 280, 20231-20238.
- Marengo E et al. (2004). Study of proteomic changes associated with healthy and tumoral murine samples in neuroblastoma by principal component analysis and classification methods. *Clinica Chimica Acta* 345, 55-67.
- Marques JM et al. (2006). *Saccharomyces cerevisiae* Hog1 Protein Phosphorylation upon Exposure to Bacterial Endotoxin. *J Biol Chem* 281, 24687-24694.
- Martin CI and Johnston IA (2005). The molecular regulation of exercised-induced muscle fibre hypertrophy in the common carp: Expression of MyoD, PCNA and components of the calcineurin-signalling pathway. *Comp Biochem Physiol Part B: Biochem Mol Biol* 142, 324-334.
- Martin CI and Johnston IA (2005). The role of myostatin and the calcineurin-signalling pathway in regulating muscle mass in response to exercise training in the rainbow trout *Oncorhynchus mykiss* Walbaum. *J Exp Biol* 208, 2083-2090.
- Masliah E et al. (2005). Effects of  $\alpha$ -Synuclein Immunization in a Mouse Model of Parkinson's Disease. *Neuron* 46, 857-868.
- Matheny CJ et al. (2004). Effect Of Prototypical Inducing Agents On P-Glycoprotein and CYP3A Expression In Mouse Tissues. *Drug Metab Dispos* 32, 1008-1014.
- Matsui H et al. (2005). Azithromycin Inhibits the Formation of Flagellar Filaments Without Suppressing Flagellin Synthesis in *Salmonella enterica* Serovar Typhimurium. *Antimicrob Agents Chemother* 49, 3396-3403.
- Matsumoto N et al. (2005). Altered cathepsin D metabolism in PHEX antisense human osteoblast cells. *Biochem Biophys Res Commun* 332, 248-253.
- Matsushima R et al. (2006). Human airway trypsin-like protease stimulates human bronchial fibroblast proliferation in a protease-activated receptor-2-dependent pathway. *Am J Physiol Lung Cell Mol Physiol* 290, L385-L395.
- McBeath R et al. (2004). Cell Shape, Cytoskeletal Tension, and RhoA Regulate Stem Cell Lineage Commitment. *Developmental Cell* 6, 483-495.
- McCarrey J (2007). Meiotic sex chromosome inactivation in the marsupial *Monodelphis domestica*. *Genesis* 45, 696-708.
- McGargill MA et al. (2005). Active Ca $^{2+}$ /Calmodulin-Dependent Protein Kinase I $\gamma$ B Impairs Positive Selection of T Cells by Modulating TCR Signaling. *J Immunol* 175, 656-664.
- Melendez L (2007). CSF proteomic fingerprints for HIV-associated cognitive impairment. *J Neuroimmunol* 192, 157-170.
- Meloni F (2007). Bronchoalveolar lavage fluid proteome in bronchiolitis obliterans syndrome: Possible role for surfactant protein A in disease onset. *J Heart Lung Transplant* 26, 1135-1143.
- Mhaidat NM et al. (2007). Docetaxel-induced apoptosis in melanoma cells is dependent on activation of caspase-2. *Mol Cancer Ther* 6, 752-761.
- Mhaidat NM et al. (2007). Docetaxel-Induced Apoptosis of Human Melanoma Is Mediated by Activation of c-Jun NH2-Terminal Kinase and Inhibited by the Mitogen-Activated Protein Kinase Extracellular Signal-Regulated Kinase 1/2 Pathway. *Clin Cancer Res* 13, 1308-1314.

- Miao J et al. (2005). Reducing Cerebral Microvascular Amyloid- $\beta$  Protein Deposition Diminishes Regional Neuroinflammation in Vasculotropic Mutant Amyloid Precursor Protein Transgenic Mice. *J Neurosci* 25, 6271-6277.
- Miglietta A et al. (2006). Conjugated linoleic acid induces apoptosis in MDA-MB-231 breast cancer cells through ERK/MAPK signalling and mitochondrial pathway. *Cancer Lett* 234, 149-157.
- Mills KV et al. (2004). Protein Splicing of a *Pyrococcus abyssi* Intein with a C-terminal Glutamine. *J Biol Chem* 279, 20685-20691.
- Milner R et al. (2007). Fibronectin- and Vitronectin-Induced Microglial Activation and Matrix Metalloproteinase-9 Expression Is Mediated by Integrins  $\alpha 5\beta 1$  and  $\alpha v\beta 5$ . *J Immunol* 178, 8158-8167.
- Milton Lima J et al. (2007). A RAPD marker identified a susceptible specific locus for gall midge resistance gene in rice cultivar ARC5984. *Crop Protection* 26, 1431-1435.
- Miyajima M et al. (2004). Effect of C-type natriuretic peptide (CNP) on water channel aquaporin-4 (AQP4) expression in cultured astrocytes. *Mol Brain Res* 122, 109-115.
- Molero JC et al. (2006). Casitas b-Lineage Lymphoma-Deficient Mice Are Protected Against High-Fat Diet-Induced Obesity and Insulin Resistance. *Diabetes* 55, 708-715.
- Molero JC et al. (2006). Genetic Ablation of the c-Cbl Ubiquitin Ligase Domain Results in Increased Energy Expenditure and Improved Insulin Action. *Diabetes* 55, 3411-3417.
- Moretti FA et al. (2007). A Major Fraction of Fibronectin Present in the Extracellular Matrix of Tissues Is Plasma-derived. *J Biol Chem* 282, 28057-28062.
- Mouécoucou J et al. (2007). Polysaccharides reduce in vitro IgG/IgE-binding of [beta]-lactoglobulin after hydrolysis. *Food Chemistry* 104, 1242-1249.
- Mulero MC et al. (2007). RCAN3, a novel calcineurin inhibitor that down-regulates NFAT-dependent cytokine gene expression. *Biochimica Biophys Acta (BBA) - Mol Cell Research* 1773, 330-341.
- Muzio G et al. (2006). Arachidonic acid suppresses growth of human lung tumor A549 cells through down-regulation of ALDH3A1 expression. *Free Rad Biol Med* 40, 1929-1938.
- Mwale F et al. (2006). The effect of glow discharge plasma surface modification of polymers on the osteogenic differentiation of committed human mesenchymal stem cells. *Biomaterials* 27, 2258-2264.
- Nagata S (2005). Isolation, characterization, and extra-embryonic secretion of the *Xenopus laevis* embryonic epidermal lectin, XEEL. *Glycobiology* 15, 281-290.
- Nakatsu Y et al. (2006). Translational Inhibition and Increased Interferon Induction in Cells Infected with C Protein-Deficient Measles Virus. *J Virol* 80, 11861-11867.
- Navarro J (2007). Effect of sublethal concentrations of copper sulphate on seabream *Sparus aurata* fingerlings. *Aquatic Living Resources* 20, 263-270.
- Nelson CM et al. (2004). Vascular Endothelial-Cadherin Regulates Cytoskeletal Tension, Cell Spreading, and Focal Adhesions by Stimulating RhoA. *Mol Biol Cell* 15, 2943-2953.
- Nezasa Ki et al. (2006). Altered Hepatobiliary Disposition Of 5 (And 6)-Carboxy-2',7'-Dichlorofluorescein In *Abcg2* (*Bcrp1*) And *Abcc2* (*Mrp2*) Knockout Mice. *Drug Metab Dispos* 34, 718-723.
- Nguyen HV et al. (2006). SUMOylation Attenuates Sensitivity Toward Hypoxia- or Desferroxamine-Induced Injury by Modulating Adaptive Responses in Salivary Epithelial Cells. *Am J Pathol* 168, 1452-1463.
- Nikolov EN and Ivanova-Nikolova TT (2007). Dynamic Integration of  $\alpha$ -Adrenergic and Cholinergic Signals in the Atria: Role Of G Protein-Regulated Inwardly Rectifying K<sup>+</sup> Channels. *J Biol Chem* 282, 28669-28682.
- Nimonkar AV and Boehmer PE (2003). The Herpes Simplex Virus Type-1 Single-strand DNA-binding Protein (ICP8) Promotes Strand Invasion. *J Biol Chem* 278, 9678-9682.
- Niture SK et al. (2007). Increased expression of the MGMT repair protein mediated by cysteine prodrugs and chemopreventative natural products in human lymphocytes and tumor cell lines. *Carcinogenesis* 28, 378-389.
- Norman RS et al. (2004). Effect of Pyocyanin on a Crude-Oil-Degrading Microbial Community. *Appl Envir Microbiol* 70, 4004-4011.
- Nyasae L et al. (2007). Dynamics of endogenous ATP7A (Menkes protein) in intestinal epithelial cells: copper-dependent redistribution between two intracellular sites. *Am J Physiol Gastrointest Liver Physiol* 292, G1181-G1194.
- Oatley JM et al. (2007). Glial Cell Line-derived Neurotrophic Factor Regulation of Genes Essential for Self-renewal of Mouse Spermatogonial Stem Cells Is Dependent on Src Family Kinase Signaling. *J Biol Chem* 282, 25842-25851.
- Oberdorster E et al. (2006). Ecotoxicology of carbon-based engineered nanoparticles: Effects of fullerene (C60) on aquatic organisms. *Carbon* 44, 1112-1120.
- Octaviani E et al. (2006). Enlazin, a Natural Fusion of Two Classes of Canonical Cytoskeletal Proteins, Contributes to Cytokinesis Dynamics. *Mol Biol Cell* 17, 5275-5286.
- Oh W et al. (2006). Jab1 Mediates Cytoplasmic Localization and Degradation of West Nile Virus Capsid Protein. *J Biol Chem* 281, 30166-30174.
- Ohnishi M et al. (2006). Integrin-linked kinase acts as a pro-survival factor against high glucose-associated osmotic stress in human mesangial cells. *Nephrol Dial Transplant* 21, 1786-1793.
- Ohno S et al. (2004). Dissection of measles virus V protein in relation to its ability to block alpha/beta interferon signal transduction. *J Gen Virol* 85, 2991-2999.
- Oliveira PJ and Wallace KB (2006). Depletion of adenine nucleotide translocator protein in heart mitochondria from doxorubicin-treated rats—Relevance for mitochondrial dysfunction. *Toxicology* 220, 160-168.
- Olson GM et al. (2007). Role of Protein O-Mannosyltransferase Pmt4 in the Morphogenesis and Virulence of *Cryptococcus neoformans*. *Eukaryot Cell* 6, 222-234.
- Overbo K and Myrnes B (2006). Deoxyribonuclease II from the Icelandic scallop (*Chlamys islandica*): Isolation and partial characterization. *Comp Biochem Physiol Part B: Biochem Mol Biol* 143, 315-318.
- Ozaslan D et al. (2003). Glycosyl Modification Facilitates Homo- and Hetero-oligomerization of the Serotonin Transporter: A Specific Role for Sialic Acid Residues. *J Biol Chem* 278, 43991-44000.
- Pandey MK et al. (2006). Induction of hepatic cytochrome P450 isozymes, benzo(a)pyrene metabolism and DNA binding following exposure to polycyclic aromatic hydrocarbon residues generated during repeated fish fried oil in rats. *Toxicol Appl Pharmacol* 213, 126-134.
- Papparella I et al. (2007). Vitamin C prevents zidovudine-induced NAD(P)H oxidase activation and hypertension in the rat. *Cardiovasc Res* 73, 432-438.
- Parmar D (2007). Overexpression of cerebral and hepatic cytochrome P450s alters behavioral activity of rat offspring following prenatal exposure to lindane. *Toxicol Appl Pharmacol* 225, 278-292.
- Patel MR et al. (2006). Effects of benzo(a)pyrene exposure on killifish (*Fundulus heteroclitus*) aromatase activities and mRNA. *Aquatic Toxicol* 77, 267-278.
- Pelley JL et al. (2007). Discovery and characterization of a novel splice variant of the GM-CSF receptor [alpha] subunit. *Exp Hematol* 35, 1483-1494.
- Pendleton AR and Machamer CE (2005). Infectious Bronchitis Virus 3a Protein Localizes to a Novel Domain of the Smooth Endoplasmic Reticulum. *J Virol* 79, 6142-6151.
- Petit A et al. (2004). Effect of cobalt and chromium ions on bcl-2, bax, caspase-3, and caspase-8 expression in human U937 macrophages. *Biomaterials* 25, 2013-2018.
- Petit A et al. (2005). Induction of protein oxidation by cobalt and chromium ions in human U937 macrophages. *Biomaterials* 26, 4416-4422.
- Petruzzella V et al. (2007). The NDUFB11 gene is not a modifier in Leber hereditary optic neuropathy. *Biochem Biophys Res Commun* 355, 181-187.
- Pfluger K and de Lorenzo V (2007). Growth-dependent Phosphorylation of the PtsN (EII<sub>Ntr</sub>) Protein of *Pseudomonas putida*. *J Biol Chem* 282, 18206-18211.
- Pham PC et al. (2004). Hypertonicity increases CLC-5 expression in mouse medullary thick ascending limb cells. *Am J Physiol Renal Physiol* 287, F747-F752.
- Phillip M (2007). Leptin regulates chondrogenic differentiation in ATDC5 cell-line through JAK/STAT and MAPK pathways. *ENDOCRINE* 32, 235-244.



- Piatek R et al. (2005). Molecular Aspects of Biogenesis of Escherichia coli Dr Fimbriae: Characterization of DraB-DraE Complexes. *Infect Immun* 73, 135-145.
- Picchi A et al. (2006). Tumor Necrosis Factor- $\alpha$  Induces Endothelial Dysfunction in the Prediabetic Metabolic Syndrome. *Circ Res* 99, 69-77.
- Piccoli C et al. (2005). Characterization of Mitochondrial and Extra-mitochondrial Oxygen Consuming Reactions in Human Hematopoietic Stem Cells: Novel Evidence of the Occurrence of NAD(P)H Oxidase Activity. *J Biol Chem* 280, 26467-26476.
- Piccoli C et al. (2007). Bone-marrow derived hematopoietic stem/progenitor cells express multiple isoforms of NADPH oxidase and produce constitutively reactive oxygen species. *Biochem Biophys Res Commun* 353, 965-972.
- Piccoli C et al. (2007). The hypoxia-inducible factor is stabilized in circulating hematopoietic stem cells under normoxic conditions. *FEBS Lett* 581, 3111-3119.
- Pindado J et al. (2007). TLR3-Dependent Induction of Nitric Oxide Synthase in RAW 264.7 Macrophage-Like Cells via a Cytosolic Phospholipase A2/Cyclooxygenase-2 Pathway. *J Immunol* 179, 4821-4828.
- Pinheiro PS et al. (2005). Presynaptic kainate receptors are localized close to release sites in rat hippocampal synapses. *Neurochem Int* 47, 309-316.
- Pinheiro PS et al. (2007). GluR7 is an essential subunit of presynaptic kainate autoreceptors at hippocampal mossy fiber synapses. *Proc Natl Acad Sci USA* 104, 12181-12186.
- Pinto-de-Oliveira A et al. (2005). Expression of HPV16 E6 oncoprotein increases resistance to several stress conditions in *Saccharomyces cerevisiae*. *FEMS Yeast Research* 5, 777-787.
- Pirone DM et al. (2006). An inhibitory role for FAK in regulating proliferation: a link between limited adhesion and RhoA-ROCK signaling. *J Cell Biol* 174, 277-288.
- Pitcovski J et al. (2005). A subunit vaccine against hemorrhagic enteritis adenovirus. *Vaccine* 23, 4697-4702.
- Place SP et al. (2004). Constitutive roles for inducible genes: evidence for the alteration in expression of the inducible hsp70 gene in Antarctic notothenioid fishes. *Am J Physiol Regulatory Integrative Comp Physiol* 287, R429-R436.
- Plante C (2007). Importance of the sedimentary matrix for anaerobic oil degradation by marine bacterial assemblages. *Bioremediation J* 11, 155-163.
- Plotkin LI et al. (2005). Bisphosphonates and Estrogens Inhibit Osteocyte Apoptosis via Distinct Molecular Mechanisms Downstream of Extracellular Signal-regulated Kinase Activation. *J Biol Chem* 280, 7317-7325.
- Plotkin LI et al. (2007). Glucocorticoids Induce Osteocyte Apoptosis by Blocking Focal Adhesion Kinase-mediated Survival: Evidence for Inside-Out Signaling Leading to Anoikis. *J Biol Chem* 282, 24120-24130.
- Plummer LJ et al. (2006). Mutational Analysis of the Ras Converting Enzyme Reveals a Requirement for Glutamate and Histidine Residues. *J Biol Chem* 281, 4596-4605.
- Polverino A et al. (2006). AMG 706, an Oral, Multikinase Inhibitor that Selectively Targets Vascular Endothelial Growth Factor, Platelet-Derived Growth Factor, and Kit Receptors, Potently Inhibits Angiogenesis and Induces Regression in Tumor Xenografts. *Cancer Res* 66, 8715-8721.
- Polymenidou M et al. (2005). Coexistence of multiple PrP<sup>Sc</sup> types in individuals with Creutzfeldt-Jakob disease. *Lancet Neurol* 4, 805-814.
- Pons S (2007). Bone morphogenetic protein 2 opposes Shh-mediated proliferation in cerebellar granule cells through a TIEG-1-based regulation of Nmyc. *J Biol Chem* 282, 37170-37180.
- Porro F et al. (2004). The erythrocyte skeletons of [beta]-adducin deficient mice have altered levels of tropomyosin, tropomodulin and EcapZ. *FEBS Lett* 576, 36-40.
- Pozzi Mucelli S et al. (2006). Proteomic studies on the white matter of human brain. *J Chromatogr B* 833, 80-90.
- Prabu SK et al. (2006). Protein Kinase A-mediated Phosphorylation Modulates Cytochrome c Oxidase Function and Augments Hypoxia and Myocardial Ischemia-related Injury. *J Biol Chem* 281, 2061-2070.
- Prost S (2007). Deficiency of G1 regulators P53, P21(Cip1) and/or pRb decreases hepatocyte sensitivity to TGF- $\beta$ ; cell cycle arrest. *BMC Cancer* 7, U1-U10.
- Provance DW, Jr. et al. (2004). From the Cover: Chemical-genetic inhibition of a sensitized mutant myosin Vb demonstrates a role in peripheral-pericentriolar membrane traffic. *Proc Natl Acad Sci USA* 101, 1868-1873.
- Puppo M et al. (2004). Induction of Apoptosis by Flavopiridol in Human Neuroblastoma Cells Is Enhanced under Hypoxia and Associated With N-myc Proto-oncogene Down-Regulation. *Clin Cancer Res* 10, 8704-8719.
- Puppo M et al. (2007). Hypoxia inhibits Moloney murine leukemia virus expression in activated macrophages. *J Leukoc Biol* 81, 528-538.
- Qian Y et al. (2003). Cloning, Characterization, and Functional Expression of the Klebsiella oxytoca XyloDEXTRIN Utilization Operon (xynTB) in Escherichia coli. *Appl Environ Microbiol* 69, 5957-5967.
- Qiu L et al. (2005). Differentially Expressed Proteins in the Pancreas of Diet-induced Diabetic Mice. *Mol Cell Proteomics* 4, 1311-1318.
- Qu J et al. (2007). Cardiac-specific haploinsufficiency of  $\beta$ -catenin attenuates cardiac hypertrophy but enhances fetal gene expression in response to aortic constriction. *J Mol Cell Cardiol* 43, 319-326.
- Rachinsky A (2007). Differential protein expression in ovaries of uninfected and Babesia-infected southern cattle ticks, Rhipicephalus (Boophilus) microplus. *Insect Biochem Mol Biol* 37, 1291-1308.
- Radovanovic I et al. (2005). Truncated Prion Protein and Doppel Are Myelinotoxic in the Absence of Oligodendrocytic PrPC. *J Neurosci* 25, 4879-4888.
- Rae C et al. (2007). Elevated NF- $\kappa$ B responses and FLIP levels in leukemic but not normal lymphocytes: reduction by salicylate allows TNF-induced apoptosis. *Proc Natl Acad Sci USA* 104, 12790-12795.
- Ramos A et al. (2007). Hypo-Osmolar Stress Induces p75NTR Expression by Activating Sp1-Dependent Transcription. *J Neurosci* 27, 1498-1506.
- Raone A et al. (2007). Hypothalamus-pituitary-adrenal modifications consequent to chronic stress exposure in an experimental model of depression in rats. *Neuroscience* 146, 1734-1742.
- Rebola N et al. (2005). Different synaptic and subsynaptic localization of adenosine A2A receptors in the hippocampus and striatum of the rat. *Neuroscience* 132, 893-903.
- Reiland J et al. (2004). Heparanase Degrades Syndecan-1 and Perlecan Heparan Sulfate: Functional Implications for Tumor Cell Invasion. *J Biol Chem* 279, 8047-8055.
- Reineri S et al. (2007). Membrane lipid rafts coordinate estrogen-dependent signaling in human platelets. *Biochimica Biophys Acta (BBA) - Mol Cell Research* 1773, 273-278.
- REN QUN et al. (2003). Comparative DNA Microarray Analysis of Host Cell Transcriptional Responses to Infection by Coxiella burnetii or Chlamydia trachomatis. *Ann N Y Acad Sci* 990, 701-713.
- Resende R et al. (2007). Susceptibility of hippocampal neurons to A $\beta$  peptide toxicity is associated with perturbation of Ca<sup>2+</sup> homeostasis. *Brain Res* 1143, 11-21.
- Reuter CJ and Maupin-Furlow JA (2004). Analysis of Proteasome-Dependent Proteolysis in Haloferax volcanii Cells, Using Short-Lived Green Fluorescent Proteins. *Appl Environ Microbiol* 70, 7530-7538.
- Richardson DK et al. (2005). Lipid Infusion Decreases the Expression of Nuclear Encoded Mitochondrial Genes and Increases the Expression of Extracellular Matrix Genes in Human Skeletal Muscle. *J Biol Chem* 280, 10290-10297.
- Richer MJ et al. (2004). Serpin Mechanism of Hepatitis C Virus Nonstructural 3 (NS3) Protease Inhibition: Induced Fit as a Mechanism For Narrow Specificity. *J Biol Chem* 279, 10222-10227.
- Richer MJ et al. (2004). The Spn4 gene of Drosophila encodes a potent furin-directed secretory pathway serpin. *Proc Natl Acad Sci USA* 101, 10560-10565.
- Richman GP et al. (2005). Quantitatively distinct requirements for signaling-competent cell spreading on engineered versus natural adhesion ligands. *J Controlled Release* 101, 3-12.
- Ripaud L et al. (2004). The [URE3] Yeast Prion Results from Protein Aggregates That Differ from Amyloid Filaments Formed in Vitro. *J Biol Chem* 279, 50962-50968.
- Robles LJ et al. (2005). Tubulin Mrna and Protein Expression are Cyclic in Retinas Exposed to Light. *Invest Ophthalmol Vis Sci* 46, 2990.

- Rockenstein E et al. (2007). Neuroprotective Effects of Regulators of the Glycogen Synthase Kinase-3 $\beta$  Signaling Pathway in a Transgenic Model of Alzheimer's Disease Are Associated with Reduced Amyloid Precursor Protein Phosphorylation. *J Neurosci* 27, 1981-1991.
- Rosca MG et al. (2005). Glycation of mitochondrial proteins from diabetic rat kidney is associated with excess superoxide formation. *Am J Physiol Renal Physiol* 289, F420-F430.
- Rossi A (2007). A quantitative and qualitative method for direct 2-DE analysis of murine cartilage. *Proteomics* 7, 4003-4007.
- Rossi A et al. (2006). Targeting the Heat Shock Factor 1 by RNA Interference: A Potent Tool to Enhance Hyperthermochemotherapy Efficacy in Cervical Cancer. *Cancer Res* 66, 7678-7685.
- Rossi D et al. (2005). Defective Tumor Necrosis Factor- $\alpha$ -dependent Control of Astrocyte Glutamate Release in a Transgenic Mouse Model of Alzheimer Disease. *J Biol Chem* 280, 42088-42096.
- Rowell C et al. (2005). Chemoprevention of Breast Cancer, Proteomic Discovery of Genistein Action in the Rat Mammary Gland. *J Nutr* 135, 2953S-2959.
- Rucevic M et al. (2006). Use of short monolithic columns for isolation of low abundance membrane proteins. *J Chromatogr A* 1123, 199-204.
- Ruiz-Manzano A et al. (2005). Levels and Activity of the Pseudomonas putida Global Regulatory Protein Crc Vary According to Growth Conditions. *J Bacteriol* 187, 3678-3686.
- Ryman KD et al. (2005). Sindbis Virus Translation Is Inhibited by a PKR/RNase L-Independent Effector Induced by Alpha/Beta Interferon Priming of Dendritic Cells. *J Virol* 79, 1487-1499.
- Sachadyn P et al. (2007). A bifunctional chimeric protein consisting of MutS and beta-galactosidase. *J Biotech* 127, 229-234.
- Saito Y et al. (2004). Synthesis and properties of novel base-discriminating fluorescent (BDF) nucleosides: a highly polarity-sensitive fluorophore for SNP typing. *Tetrahedron Lett* 45, 7827-7831.
- Sakamoto A (2007). The RNAi-mediated silencing of xanthine dehydrogenase impairs growth and fertility and accelerates leaf senescence in transgenic Arabidopsis plants. *Plant Cell Physiol* 48, 1484-1495.
- Sakamoto A et al. (2004). Three distinct Arabidopsis hemoglobins exhibit peroxidase-like activity and differentially mediate nitrite-dependent protein nitration. *FEBS Lett* 572, 27-32.
- Sanchez-Martin CC et al. (2007). Cholesterol Starvation Induces Differentiation of Human Leukemia HL-60 Cells. *Cancer Res* 67, 3379-3386.
- Sancho R et al. (2004). Imperatorin Inhibits HIV-1 Replication through an Sp1-dependent Pathway. *J Biol Chem* 279, 37349-37359.
- Sancho R et al. (2005). Mechanisms of HIV-1 Inhibition by the Lipid Mediator N-Arachidonoyldopamine. *J Immunol* 175, 3990-3999.
- Santiago AR et al. (2007). High glucose induces caspase-independent cell death in retinal neural cells. *Neurobiol Dis* 25, 464-472.
- Saoncella S et al. (2004). Syndecan-4 Regulates ATF-2 Transcriptional Activity in a Rac1-dependent Manner. *J Biol Chem* 279, 47172-47176.
- Sardanelli AM et al. (2006). Occurrence of A-kinase anchor protein and associated cAMP-dependent protein kinase in the inner compartment of mammalian mitochondria. *FEBS Lett* 580, 5690-5696.
- Sartori M et al. (2004). Effects of angiotensin II and insulin on ERK1/2 activation in fibroblasts from hypertensive patients. *Ame J Hypertension* 17, 604-610.
- Sathaye A et al. (2006). Electrical pacing counteracts intrinsic shortening of action potential duration of neonatal rat ventricular cells in culture. *J Mol Cell Cardiol* 41, 633-641.
- Sato Y et al. (2007). Stable integration and conditional expression of electroporated transgenes in chicken embryos. *Dev Biol* 305, 616-624.
- Sawhney RS et al. (2006). Integrin  $\alpha$ 2-mediated ERK and Calpain Activation Play a Critical Role in Cell Adhesion and Motility via Focal Adhesion Kinase Signaling: Identification of a Novel Signaling Pathway. *J Biol Chem* 281, 8497-8510.
- Sbodio JI and Machamer CE (2007). Identification of a Redox-sensitive Cysteine in GCP60 That Regulates Its Interaction with Golgin-160. *J Biol Chem* 282, 29874-29881.
- Sbodio JI et al. (2006). GCP60 Preferentially Interacts with a Caspase-generated Golgin-160 Fragment. *J Biol Chem* 281, 27924-27931.
- Scacco S et al. (2003). Pathological Mutations of the Human NDUFS4 Gene of the 18-kDa (A/QDQ) Subunit of Complex I Affect the Expression of the Protein and the Assembly and Function of the Complex. *J Biol Chem* 278, 44161-44167.
- Schauble S et al. (2007). Identification of ChChd3 as a Novel Substrate of the cAMP-dependent Protein Kinase (PKA) Using an Analog-sensitive Catalytic Subunit. *J Biol Chem* 282, 14952-14959.
- Schild A et al. (2006). Altered levels of PP2A regulatory B/PR55 isoforms indicate role in neuronal differentiation. *International J Dev Neurosci* 24, 437-443.
- Schwanninger R et al. (2007). Lack of Noggin Expression by Cancer Cells Is a Determinant of the Osteoblast Response in Bone Metastases. *Am J Pathol* 170, 160-175.
- See J et al. (2004). Oligodendrocyte maturation is inhibited by bone morphogenetic protein. *Mol Cell Neurosci* 26, 481-492.
- Semba S and Huebner K (2006). Protein Expression Profiling Identifies Cyclophilin A as a Molecular Target in Fhit-Mediated Tumor Suppression. *Mol Cancer Res* 4, 529-538.
- Seol GH et al. (2007). Neuromodulators Control the Polarity of Spike-Timing-Dependent Synaptic Plasticity. *Neuron* 55, 919-929.
- Serviddio G et al. (2005). Brief hypoxia before normoxic reperfusion (postconditioning) protects the heart against ischemia-reperfusion injury by preventing mitochondrial peroxide production and glutathione depletion. *FASEB J* 19, 354-361.
- Shadel G (2007). Convergence of multiple signaling pathways is required to coordinately up-regulate mtDNA and mitochondrial biogenesis during T cell activation. *Mitochondrion* 7, 374-385.
- Shearer J et al. (2005). Partial Gene Deletion of Heart-Type Fatty Acid-Binding Protein Limits the Severity of Dietary-Induced Insulin Resistance. *Diabetes* 54, 3133-3139.
- Shen Z (2007). Bacillus thuringiensis Vip3 mutant proteins: Insecticidal activity and trypsin sensitivity. *Biocontrol Sci Technol* 17, 699-708.
- Shi L et al. (2007). Caloric restriction eliminates the aging-related decline in NMDA and AMPA receptor subunits in the rat hippocampus and induces homeostasis. *Exp Neurol* 206, 70-79.
- Shutman M et al. (2006). Cell Adhesion Molecule L1 Disrupts E-Cadherin-Containing Adherens Junctions and Increases Scattering and Motility of MCF7 Breast Carcinoma Cells. *Cancer Res* 66, 11370-11380.
- Shukla Y (2007). In vitro and in vivo modulation of testosterone mediated alterations in apoptosis related proteins by [6]-gingerol. *Mol Nutri Food Res* 51, 1492-1502.
- Shukla Y (2007). Theaflavins induce G2/M arrest by modulating expression of p21(waf1/cip1), cdc25C and cyclin B in human prostate carcinoma PC-3 cells. *Life Sci* 81, 1323-1331.
- Shults CW et al. (2005). Neurological and Neurodegenerative Alterations in a Transgenic Mouse Model Expressing Human  $\alpha$ -Synuclein under Oligodendrocyte Promoter: Implications for Multiple System Atrophy. *J Neurosci* 25, 10689-10699.
- Si Y et al. (2007). Effects of forced uncoupling protein 1 expression in 3T3-L1 cells on mitochondrial function and lipid metabolism. *J Lipid Res* 48, 826-836.
- Siciliano V et al. (2007). Transcriptome Analysis of Arbuscular Mycorrhizal Roots during Development of the Prepenetration Apparatus. *Plant Physiol* 144, 1455-1466.
- Sigdel TK et al. (2004). Fractionation of Soluble Proteins in Escherichia coli Using DEAE-, SP-, and Phenyl Sepharose Chromatographies. *J Biomol Tech* 15, 199-207.
- Silva CG et al. (2007). Blockade of adenosine A2A receptors prevents staurosporine-induced apoptosis of rat hippocampal neurons. *Neurobiol Dis* 27, 182-189.
- Silva VS et al. (2005). Effect of Chronic Exposure to Aluminium on Isoform Expression and Activity of Rat (Na+/K+) ATPase. *Toxicol Sci* 88, 485-494.
- Simon R and Samuel CE (2007). Activation of NF- $\kappa$ B-dependent gene expression by Salmonella flagellins FliC and FliB. *Biochem Biophys Res Commun* 355, 280-285.

- Slominski A et al. (2005). CRH stimulates POMC activity and corticosterone production in dermal fibroblasts. *J Neuroimmunol* 162, 97-102.
- Smallwood ND et al. (2005). Involvement of NH2 terminus of PKC- $\delta$  in binding to F-actin during activation of Calu-3 airway epithelial NKCC1. *Am J Physiol Cell Physiol* 288, C906-C912.
- Stanislawska-Sachadyn A et al. (2005). Preliminary studies on DNA retardation by MutS applied to the detection of point mutations in clinical samples. *Mutation Research/Fundamental and Molecular Mechanisms of Mutagenesis* 570, 97-103.
- Stanislawska-Sachadyn A et al. (2006). The construction of bifunctional fusion proteins consisting of MutS and GFP. *J Biotech* 121, 134-143.
- Stefani G et al. (2004). Fragile X Mental Retardation Protein Is Associated with Translating Polyribosomes in Neuronal Cells. *J Neurosci* 24, 7272-7276.
- Steinhauer D (2007). Analysis of residues near the fusion peptide in the influenzahemagglutinin structure for roles in triggering membrane fusion. *Virology* 370, 403-414.
- Stewart D et al. (2004). Ubiquitination and proteasome degradation of the E6 proteins of human papillomavirus types 11 and 18. *J Gen Virol* 85, 1419-1426.
- Stocum ET and Plante CJ (2006). The effect of artificial defaunation on bacterial assemblages of intertidal sediments. *J Exp Mar Biol Ecol* 337, 147-158.
- Su D (2007). The aged thymus shows normal recruitment of lymphohematopoietic progenitors but has defects in thymic epithelial cells. *Int Immunol* 19, 1201-1211.
- Suchankova G et al. (2005). Dietary polyunsaturated fatty acids enhance hepatic AMP-activated protein kinase activity in rats. *Biochem Biophys Res Commun* 326, 851-858.
- Sunol C (2007). Long-term exposure to dieldrin reduces  $\gamma$ -aminobutyric acid type A and N-methyl-D-aspartate receptor function in primary cultures of mouse cerebellar granule cells. *J Neurosci Res* 85, 3687-3695.
- Sutton RB et al. (2005). Crystal Structure of Cone Arrestin at 2.3 Å: Evolution of Receptor Specificity. *J Mol Biol* 354, 1069-1080.
- Suzuki K et al. (2006). Identification of a novel regulatory protein (CsrD) that targets the global regulatory RNAs CsrB and CsrC for degradation by RNase E. *Genes and Dev* 20, 2605-2617.
- Szweda P (2007). Molecular typing of *Staphylococcus aureus* isolated from cows with mastitis in the east of Poland on the basis of polymorphism of genes coding protein A and coagulase. *Polish J Vet Sci* 10, 199-205.
- Tahara M et al. (2007). Altered Interaction of the Matrix Protein with the Cytoplasmic Tail of Hemagglutinin Modulates Measles Virus Growth by Affecting Virus Assembly and Cell-Cell Fusion. *J Virol* 81, 6827-6836.
- Tain YL et al. (2007). Vitamin E reduces glomerulosclerosis, restores renal neuronal NOS, and suppresses oxidative stress in the 5/6 nephrectomized rat. *Am J Physiol Renal Physiol* 292, F1404-F1410.
- Takeda M et al. (2005). Long Untranslated Regions of the Measles Virus M and F Genes Control Virus Replication and Cytopathogenicity. *J Virol* 79, 14346-14354.
- Talarek N et al. (2005). The [URE3] Prion Is Not Conserved Among *Saccharomyces* Species. *Genetics* 171, 23-34.
- Talarico LA et al. (2005). Construction and expression of an ethanol production operon in Gram-positive bacteria. *Microbiology* 151, 4023-4031.
- Tallman MN et al. (2007). The Contribution of Intestinal UDP-Glucuronosyltransferases in Modulating 7-Ethyl-10-hydroxy-camptothecin (SN-38)-Induced Gastrointestinal Toxicity in Rats. *J Pharmacol Exp Ther* 320, 29-37.
- Tang ES and Martin A (2006). NHEJ-deficient DT40 cells have increased levels of immunoglobulin gene conversion: evidence for a double strand break intermediate. *Nucleic Acids Res* 34, 6345-6351.
- Tang X et al. (2005). LPS induces the interaction of a transcription factor, LPS-induced TNF- $\alpha$  factor, and STAT6(B) with effects on multiple cytokines. *Proc Natl Acad Sci USA* 102, 5132-5137.
- Tang X et al. (2006). LPS-induced TNF- $\alpha$  factor (LITAF)-deficient mice express reduced LPS-induced cytokine: Evidence for LITAF-dependent LPS signaling pathways. *Proc Natl Acad Sci USA* 103, 13777-13782.
- Tang YD et al. (2005). Low Thyroid Function Leads to Cardiac Atrophy With Chamber Dilatation, Impaired Myocardial Blood Flow, Loss of Arterioles, and Severe Systolic Dysfunction. *Circulation* 112, 3122-3130.
- Tao RH et al. (2006). Testicular zinc finger protein recruits histone deacetylase 2 and suppresses the transactivation function and intranuclear foci formation of agonist-bound androgen receptor competitively with TIF2. *Mol Cell Endocrinol* 247, 150-165.
- Tardito D et al. (2007). Long-term soluble A $\beta$ 1-40 activates CaM kinase II in organotypic hippocampal cultures. *Neurobiol Aging* 28, 1388-1395.
- Taru H et al. (2004). Suppression of the caspase cleavage of  $\beta$ -amyloid precursor protein by its cytoplasmic phosphorylation. *FEBS Lett* 567, 248-252.
- Taverna E et al. (2004). Role of Lipid Microdomains in P/Q-type Calcium Channel (Cav2.1) Clustering and Function in Presynaptic Membranes. *J Biol Chem* 279, 5127-5134.
- Temmink OH et al. (2007). Irinotecan-induced cytotoxicity to colon cancer cells in vitro is stimulated by pre-incubation with trifluorothymidine. *Euro J Cancer* 43, 175-183.
- Teodoro J et al. (2006). Decreased ANT content in Zucker fatty rats: Relevance for altered hepatic mitochondrial bioenergetics in steatosis. *FEBS Lett* 580, 2153-2157.
- Tesco G et al. (2007). Depletion of GGA3 Stabilizes BACE and Enhances  $\beta$ -Secretase Activity. *Neuron* 54, 721-737.
- Thomas JM et al. (2003). Sindbis Virus Vectors Designed To Express a Foreign Protein as a Cleavable Component of the Viral Structural Polyprotein. *J Virol* 77, 5598-5606.
- Thomas TA et al. (2005). Thyroid hormones induce unique and potentially beneficial changes in cardiac myocyte shape in hypertensive rats near heart failure. *Am J Physiol Heart Circ Physiol* 288, H2118-H2122.
- Ting HJ et al. (2007). Docetaxel-induced growth inhibition and apoptosis in androgen independent prostate cancer cells are enhanced by 1 $\alpha$ ,25-dihydroxyvitamin D3. *Cancer Lett* 247, 122-129.
- Tjwe SL et al. (2005). Expression of N-acetylglucosamine 6-O-sulfotransferases (GlcNAc6STs)-1 and -4 in human monocytes: GlcNAc6ST-1 is implicated in the generation of the 6-sulfo N-acetylglucosamine/Lewis x epitope on CD44 and is induced by TNF- $\alpha$ . *Glycobiology* 15, 7C-13.
- Toporowski MC et al. (2004). Two Outer Membrane Proteins Are Required for Maximal Type I Secretion of the *Caulobacter crescentus* S-Layer Protein. *J Bacteriol* 186, 8000-8009.
- Torrado M et al. (2004). Left-right asymmetric ventricular expression of CARP in the piglet heart: regional response to experimental heart failure. *Euro J Heart Failure* 6, 161-172.
- Turncliff RZ et al. (2004). Effect Of Dexamethasone Treatment On The Expression And Function Of Transport Proteins In Sandwich-Cultured Rat Hepatocytes. *Drug Metab Dispos* 32, 834-839.
- Umemiya R et al. (2007). Cloning and characterization of an autophagy-related gene, ATG12, from the three-host tick *Haemaphysalis longicornis*. *Insect Biochem Mol Biol* 37, 975-984.
- Umemiya R et al. (2007). *Haemaphysalis longicornis*: Molecular characterization of a homologue of the macrophage migration inhibitory factor from the partially fed ticks. *Exp Parasitol* 115, 135-142.
- Uziel O et al. (2007). Telomere dynamics in arteries and mononuclear cells of diabetic patients: Effect of diabetes and of glycemic control. *Exp Gerontol* 42, 971-978.
- Vaccari G et al. (2006). Identification of an allelic variant of the goat PrP gene associated with resistance to scrapie. *J Gen Virol* 87, 1395-1402.
- Van Nostrand W (2007). Induction of complement proteins in a mouse model for cerebral microvascular A $\beta$  deposition. *J Neuroinflammation* 4, U1-U8.
- Varo I et al. (2007). Effects of dichlorvos aquaculture treatments on selected biomarkers of gilthead sea bream (*Sparus aurata* L.) fingerlings. *Aquaculture* 266, 87-96.
- Vascotto F et al. (2004). Effects of intrabodies specific for rotavirus NSP5 during the virus replicative cycle. *J Gen Virol* 85, 3285-3290.

- Venkatakrishnan CD et al. (2006). Heat shock protects cardiac cells from doxorubicin-induced toxicity by activating p38 MAPK and phosphorylation of small heat shock protein 27. *Am J Physiol Heart Circ Physiol* 291, H2680-H2691.
- Vidali M et al. (2007). Detection of auto-antibodies against cytochrome P4502E1 (CYP2E1) in chronic hepatitis C. *J Hepatology* 46, 605-612.
- Vincent D et al. (2007). Proteomic analysis reveals differences between Vitis vinifera L. cv. Chardonnay and cv. Cabernet Sauvignon and their responses to water deficit and salinity. *J Exp Bot* 58, 1873-1892.
- Vollrath D (2007). Rapid and stable knockdown of an endogenous gene in retinal pigment epithelium. *Hum Gene Ther* 18, 871-880.
- Voltan S et al. (2007). Aggregating Phenotype in *Lactobacillus crispatus* Determines Intestinal Colonization and TLR2 and TLR4 Modulation in Murine Colonic Mucosa. *Clin Vaccine Immunol* 14, 1138-1148.
- Wang D et al. (2006). Systemic aspects of host-pathogen interactions in Austrian pine (*Pinus nigra*): A proteomics approach. *Physiol Mol Plant Pathol* 68, 149-157.
- Wang H et al. (2005). DNA Ligase III as a Candidate Component of Backup Pathways of Nonhomologous End Joining. *Cancer Res* 65, 4020-4030.
- Wang M (2007). Blocking tumorigenic activities of colorectal cancer cells by asplicing RON receptor variant defective in the tyrosine kinase domain. *Cancer Biol Ther* 6, 1121-1129.
- Wang P et al. (2007). Physical and Functional Interactions of SNAP-23 with Annexin A2. *Am J Respir Cell Mol Biol* 37, 467-476.
- Wang X et al. (2007). Ethanol attenuates Aldh9 mRNA expression in Japanese medaka (*Oryzias latipes*) embryogenesis. *Comp Biochem Physiol Part B: Biochem Mol Biol* 146, 357-363.
- Wang YF et al. (2007). Apoptosis Induction in Human Melanoma Cells by Inhibition of MEK Is Caspase-Independent and Mediated by the Bcl-2 Family Members PUMA, Bim, and Mcl-1. *Clin Cancer Res* 13, 4934-4942.
- Wang Z (2007). High efficient expression and bio-activity assay of recombinant antibody for citrus bacterial canker disease. *Weishengwu Xuebao* 47, 1066-1069.
- Wang Z et al. (2007). Human Mitochondrial Ribosomal Protein MRPL12 Interacts Directly with Mitochondrial RNA Polymerase to Modulate Mitochondrial Gene Expression. *J Biol Chem* 282, 12610-12618.
- Wehkamp J et al. (2005). Reduced Paneth cell  $\alpha$ -defensins in ileal Crohn's disease. *Proc Natl Acad Sci USA* 102, 18129-18134.
- Wei W et al. (2006). The LDL Receptor-Related Protein LRP6 Mediates Internalization and Lethality of Anthrax Toxin. *Cell* 124, 1141-1154.
- Weitsman GE et al. (2005). Vitamin D sensitizes breast cancer cells to the action of H<sub>2</sub>O<sub>2</sub>: Mitochondria as a convergence point in the death pathway. *Free Rad Biol Med* 39, 266-278.
- Wen VW et al. (2006). Telomere-Driven Karyotypic Complexity Concur with p16INK4a Inactivation in TP53-Competent Immortal Endothelial Cells. *Cancer Res* 66, 10691-10700.
- Wente W et al. (2006). Fibroblast Growth Factor-21 Improves Pancreatic  $\beta$ -Cell Function and Survival by Activation of Extracellular Signal-Regulated Kinase 1/2 and Akt Signaling Pathways. *Diabetes* 55, 2470-2478.
- Wente W et al. (2007). Activation of Liver X Receptors and Retinoid X Receptors Induces Growth Arrest and Apoptosis in Insulin-Secreting Cells. *Endocrinology* 148, 1843-1849.
- White NM et al. (2004). Mechanistic Similarities between Cultured Cell Models of Cystic Fibrosis and Niemann-Pick Type C. *Am J Respir Cell Mol Biol* 31, 538-543.
- White NM et al. (2007). Altered cholesterol homeostasis in cultured and in vivo models of cystic fibrosis. *Am J Physiol Lung Cell Mol Physiol* 292, L476-L486.
- Whitlock NA et al. (2005). Hsp27 Upregulation by HIF-1 Signaling Offers Protection against Retinal Ischemia in Rats. *Invest Ophthalmol Vis Sci* 46, 1092-1098.
- Williams SK et al. (2007). A novel epiphytic cyanobacterium associated with reservoirs affected by avian vacuolar myelinopathy. *Harmful Algae* 6, 343-353.
- Woo JL and Berk AJ (2007). Adenovirus Ubiquitin-Protein Ligase Stimulates Viral Late mRNA Nuclear Export. *J Virol* 81, 575-587.
- Xia Y and Krukoff TL (2004). Estrogen Induces Nitric Oxide Production via Activation of Constitutive Nitric Oxide Synthases in Human Neuroblastoma Cells. *Endocrinology* 145, 4550-4557.
- Xu F et al. (2005). Protease nexin-2/amyloid  $\beta$ -protein precursor limits cerebral thrombosis. *Proc Natl Acad Sci USA* 102, 18135-18140.
- Xu F et al. (2007). Early-onset subicular microvascular amyloid and neuroinflammation correlate with behavioral deficits in vasculotropic mutant amyloid  $\beta$ -protein precursor transgenic mice. *Neuroscience* 146, 98-107.
- Xu S and Hori RT (2004). Identification of a domain within human TAF<sub>148</sub>, a subunit of Selectivity Factor 1, that interacts with helix 2 of TBP. *Gene* 338, 177-186.
- Xu X et al. (2006). Shared Long-Range Regulatory Elements Coordinate Expression of a Gene Cluster Encoding Nicotinic Receptor Heteromeric Subtypes. *Mol Cell Biol* 26, 5636-5649.
- Xu X et al. (2007). Anti-LOX-1 Rescues Endothelial Function in Coronary Arterioles in Atherosclerotic ApoE Knockout Mice. *Arterioscler Thromb Vasc Biol* 27, 871-877.
- Yadav S et al. (2006). Regional specificity in deltamethrin induced cytochrome P450 expression in rat brain. *Toxicol Appl Pharmacol* 217, 15-24.
- Yamada O et al. (2007). Light-chain multiple myeloma in a cat. *J Vet Diagn Invest* 19, 443-447.
- Yamaguchi K et al. (2006). Epidermal growth factor-induced mobilization of a ganglioside-specific sialidase (NEU3) to membrane ruffles. *Biochem Biophys Res Commun* 346, 484-490.
- Yamamoto-Mizuma S et al. (2004). Altered properties of volume-sensitive osmolyte and anion channels (VSOACs) and membrane protein expression in cardiac and smooth muscle myocytes from *Clcn3*<sup>-/-</sup> mice. *J Physiol* 557, 439-456.
- Yamazaki-Sato H et al. (2003). The carboxyl terminal 17 amino acids within Apg7 are essential for Apg8 lipidation, but not for Apg12 conjugation. *FEBS Lett* 551, 71-77.
- Yang R and Newgard CB (2003). Hepatic Expression of a Targeting Subunit of Protein Phosphatase-1 in Streptozotocin-diabetic Rats Reverses Hyperglycemia and Hyperphagia Despite Depressed Glucokinase Expression. *J Biol Chem* 278, 23418-23425.
- Yang Y and Loscalzo J (2005). S-nitrosoprotein formation and localization in endothelial cells. *Proc Natl Acad Sci USA* 102, 117-122.
- Yao H (2007). Altered expression of the RON receptor tyrosine kinase in various epithelial cancers and its contribution to tumorigenic phenotypes in thyroid cancer cells. *J Pathol* 213, 402-411.
- Yarian CS et al. (2005). Aconitase and ATP synthase are targets of malondialdehyde modification and undergo an age-related decrease in activity in mouse heart mitochondria. *Biochem Biophys Res Commun* 330, 151-156.
- Yeh S et al. (2003). Abnormal Mammary Gland Development and Growth Retardation in Female Mice and MCF7 Breast Cancer Cells Lacking Androgen Receptor. *J Exp Med* 198, 1899-1908.
- Yeom JH and Lee K (2006). RraA rescues *Escherichia coli* cells over-producing RNase E from growth arrest by modulating the ribonucleolytic activity. *Biochem Biophys Res Commun* 345, 1372-1376.
- Yeom SY et al. (2006). Regulation of Insulin Secretion and  $\beta$ -Cell Mass by Activating Signal Cointegrator 2. *Mol Cell Biol* 26, 4553-4563.
- Ying WZ and Sanders PW (2005). Enhanced expression of EGF receptor in a model of salt-sensitive hypertension. *Am J Physiol Renal Physiol* 289, F314-F321.
- Ying WZ et al. (2007). EGF Receptor Activity Modulates Apoptosis Induced by Inhibition of the Proteasome of Vascular Smooth Muscle Cells. *J Am Soc Nephrol* 18, 131-142.
- Yu D (2007). Comparative proteomic analysis of wild (Glycine soja) and cultivated (Glycine max) soybean seeds. *Prog Inorg Biochem Biophys* 34, 1296-1302.



- Yu Z and Shah DM (2007). Curcumin down-regulates Ets-1 and Bcl-2 expression in human endometrial carcinoma HEC-1-A cells. *Gynecologic Oncology* 106, 541-548.
- Yue H et al. (2007). Effect of nitrogen application rate on content of glutenin macropolymer and high molecular weight glutenin subunits in grains of two winter wheat cultivars. *J Cereal Sci* 45, 248-256.
- Yuhas Y (2007). Rifampin inhibits prostaglandin E(2) production and arachidonic acid release in human alveolar epithelial cells. *Antimicrob Agents Chemother* 51, 4225-4230.
- Yuhas Y et al. (2006). Rifampin Augments Cytokine-Induced Nitric Oxide Production in Human Alveolar Epithelial Cells. *Antimicrob Agents Chemother* 50, 396-398.
- Yung E et al. (2004). Specificity of Interaction of IN1/hSNF5 with Retroviral Integrases and Its Functional Significance. *J Virol* 78, 2222-2231.
- Zalewska B et al. (2005). A surface-exposed DraD protein of uropathogenic *Escherichia coli* bearing Dr fimbriae may be expressed and secreted independently from DraC usher and DraE adhesin. *Microbiology* 151, 2477-2486.
- Zarembek V et al. (2005). Cytotoxicity of an Anti-cancer Lysophospholipid through Selective Modification of Lipid Raft Composition. *J Biol Chem* 280, 38047-38058.
- Zeman J (2007). Knockdown of human Oxa1l impairs the biogenesis of F(1)F(o)-ATP synthase and NADH:Ubiquinone oxidoreductase. *J Mol Biol* 374, 506-516.
- Zeng W (2007). Temporal and spatial changes of histone 3 K4 dimethylation at the FN- $\gamma$  gene during Th1 and Th2 cell differentiation. *J Immunol* 179, 6410-6415.
- Zhang P and Samuel CE (2007). Protein Kinase PKR Plays a Stimulus- and Virus-Dependent Role in Apoptotic Death and Virus Multiplication in Human Cells. *J Virol* 81, 8192-8200.
- Zhang Y et al. (2005). Adenosine-Dependent Induction of Glutathione Peroxidase 1 in Human Primary Endothelial Cells and Protection Against Oxidative Stress. *Circ Res* 96, 831-837.
- Zhang Y et al. (2007). Identification and Characterization of Interferon-Induced Proteins That Inhibit Alphavirus Replication. *J Virol* 81, 11246-11255.
- Zhang YJ et al. (2007). Inhibition of replication and transcription activator and latency-associated nuclear antigen of Kaposi's sarcoma-associated herpesvirus by morpholino oligomers. *Antiviral Res* 73, 12-23.
- Zheng G and Yang YC (2004). ZNF76, a Novel Transcriptional Repressor Targeting TATA-binding Protein, Is Modulated by Sumoylation. *J Biol Chem* 279, 42410-42421.
- Zhou J et al. (2007). Upregulation of  $\gamma$ -catenin compensates for the loss of beta-catenin in adult cardiomyocytes. *Am J Physiol Heart Circ Physiol* 292, H270-H276.
- Zhou Q and Amar S (2006). Identification of Proteins Differentially Expressed in Human Monocytes Exposed to *Porphyromonas gingivalis* and Its Purified Components by High-Throughput Immunoblotting. *Infect Immun* 74, 1204-1214.
- Zhou Q et al. (2005). Cytokine Profiling of Macrophages Exposed to *Porphyromonas gingivalis*, Its Lipopolysaccharide, or Its FimA Protein. *Infect Immun* 73, 935-943.
- Zhu J et al. (2003). Posttranscriptional control of the competence factor  $\beta$ FTZ-F1 by juvenile hormone in the mosquito *Aedes aegypti*. *Proc Natl Acad Sci USA* 100, 13338-13343.
- Zhu Qs et al. (2004). CCAAT/Enhancer-binding Protein  $\alpha$  (C/EBP $\alpha$ ) Activates Transcription of the Human Microsomal Epoxide Hydrolase Gene (EPHX1) through the Interaction with DNA-bound NF- $\kappa$ B. *J Biol Chem* 279, 29902-29910.
- Zhu Qs et al. (2004). Regulation of human microsomal epoxide hydrolase gene (EPHX1) expression by the transcription factor GATA-4. *Biochimica Biophys Acta (BBA) – Gene Structure and Expression* 1676, 251-260.
- Zhu S et al. (2006). Environmental induction of CYP1A-, CYP2M1- and CYP2K1-like proteins in tropical fish species by produced formation water on the northwest shelf of Australia. *Marine Environ Res* 62, S322-S326.
- Zianni M et al. (2006). Identification of the DNA Bases of a DNase I Footprint by the Use of Dye Primer Sequencing on an Automated Capillary DNA Analysis Instrument. *J Biomol Tech* 17, 103-113.
- Zimman A et al. (2007). Vascular Endothelial Growth Factor Receptor 2 Plays a Role in the Activation of Aortic Endothelial Cells by Oxidized Phospholipids. *Arterioscler Thromb Vasc Biol* 27, 332-338.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

---

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400  
**Canada** 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65  
**Germany** 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **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** 0508 805 500  
**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** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000

---