

Quantity One® 1-D Analysis Software Citations

- Arien KK, Abraha A, Quinones-Mateu ME, Kestens L, Vanham G, and Arts EJ, The replicative fitness of primary human immunodeficiency virus type 1 (HIV-1) group M, HIV-1 group O, and HIV-2 isolates, *J Virol* 79, 8979–8990 (2005)
- Bandiera S, Weidlich S, Harth V, Broede P, Ko Y, and Friedberg T, Proteasomal degradation of human CYP1B1: effect of the Asn453Ser polymorphism on the post-translational regulation of CYP1B1 expression, *Mol Pharmacol* 67, 435–443 (2005)
- Carbone MA, Llopart A, deAngelis M, Coyne JA, and Mackay TFC, Quantitative trait loci affecting the difference in pigmentation between *Drosophila yakuba* and *D. santomea*, *Genetics* 171, 211–225 (2005)
- Choi JY and Guengerich FP, Kinetic evidence for inefficient and error-prone bypass across bulky N₂-guanine DNA adducts by human DNA polymerase, *J Biol Chem*, M600112200, published electronically March 8 (2006)
- Dorweiler IJ, Ruone SJ, Wang H, Burry RW, and Mansky LM, Role of the human T-cell leukemia virus type 1 PTAP motif in Gag targeting and particle release, *J Virol* 80, 3634–3643 (2006)
- Ekman M, Fagher K, Wede M, Stakeberg K, and Arner A, Decreased phosphatase activity, increased Ca²⁺ sensitivity, and myosin light chain phosphorylation in urinary bladder smooth muscle of newborn mice, *J Gen Physiol* 125, 187–196 (2005)
- Enwall K, Philippot L, and Hallin S, Activity and composition of the denitrifying bacterial community respond differently to long-term fertilization, *Appl Environ Microbiol* 71, 8335–8343 (2005)
- Fiola K, Perreault JP, and Cousineau B, Gene targeting in the gram-positive bacterium *Lactococcus lactis*, using various delta ribozymes, *Appl Environ Microbiol* 72, 869–879 (2006)
- Han ES, Cooper DL, Persky NS, Sutura VA Jr, Whitaker RD, Montello ML, and Lovett ST, RecJ exonuclease: substrates, products and interaction with SSB, *Nucleic Acids Res* 34, 1084–1091 (2006)
- Jimenez J, Jang GM, Semler BL, and Waterman ML, An internal ribosome entry site mediates translation of lymphoid enhancer factor-1, *RNA* 11, 1385–1399 (2005)
- Katsiari CG, Kytтары VC, Juang YT, and Tsokos GC, Protein phosphatase 2A is a negative regulator of IL-2 production in patients with systemic lupus erythematosus, *J Clin Invest* 115, 3193–3204 (2005)
- Kinlough CL, McMahan RJ, Poland PA, Bruns JB, Harkleroad KL, Stremple RJ, Kashlan OB, Weixel KM, Weisz OA, and Hughey RP, MUC1 recycling is dependent on its palmitoylation, *J Biol Chem*, M512996200, published electronically February 28 (2006)
- Liu J, Chen Q, Huang W, Horak KM, Zheng H, Mestrl R, and Wang X, Impairment of the ubiquitin-proteasome system in desminopathy mouse hearts, *FASEB J* 2, 362–364 (2005)
- Liu Q, White LR, Clark SA, Heffner DJ, Winston BW, Tibbles LA, and Muruve DA, Akt/protein kinase B activation by adenovirus vectors contributes to NFκB-dependent CXCL10 expression, *J Virol* 79, 14507–14515 (2005)
- McCabe, CD and Innis JW, A genomic approach to the identification and characterization of HOXA13 functional binding elements, *Nucleic Acids Res* 33, 6782–6794 (2005)
- Mezosi E, Wang SH, Utsugi S, Bajnok L, Bretz JD, Gauger PG, Thompson NW, and Baker JR Jr, Induction and regulation of Fas-mediated apoptosis in human thyroid epithelial cells, *Mol Endocrinol* 19, 804–811 (2005)
- Moreau K, Dizin E, Ray H, Luquain C, Lefai E, Foufelle F, Billaud M, Lenoir GM, and Venezia ND, BRCA1 affects lipid synthesis through its interaction with acetyl-CoA carboxylase, *J Biol Chem* 281, 3172–3181 (2006)
- Rahayu YS, Walch-Liu P, Neumann G, Romheld V, von Wiren N, and Bangerth F, Root-derived cytokinins as long-distance signals for NO₃⁻-induced stimulation of leaf growth, *J Exp Bot* 56, 1143–1152 (2005)
- Rosenkilde MM, Benned-Jensen T, Andersen H, Holst PJ, Kledal TN, Luttichau HR, Larsen JK, Geisler C, Christensen JP, and Schwartz TW, Molecular pharmacological phenotyping of EBI2 — an orphan 7TM receptor with constitutive activity, *J Biol Chem*, M602245200, published electronically March 15 (2006)
- Seong IS, Ivanova E, Lee JM, Choo YS, Fossale E, Anderson M, Gusella JF, Laramie JM, Myers RH, Lesort M, and MacDonald ME, HD CAG repeat implicates a dominant property of huntingtin in mitochondrial energy metabolism, *Hum Mol Genet* 14, 2871–2880 (2005)
- Shao J, Qiao L, Janssen RC, Pagliassotti M, and Friedman JE, Chronic hyperglycemia enhances PEPCK gene expression and hepatocellular glucose production via elevated liver activating protein/liver inhibitory protein ratio, *Diabetes* 54, 976–984 (2005)
- Shay KP, Wang Z, Xing PX, McKenzie IFC, and Magnuson NS, Pim-1 kinase stability is regulated by heat shock proteins and the ubiquitin-proteasome pathway, *Mol Cancer Res* 3, 170–181 (2005)
- Sommer U, Schmid C, Sobota RM, Lehmann U, Stevenson NJ, Johnston JA, Schaper F, Heinrich PC, and Haan S, Mechanisms of SOCS3 phosphorylation upon interleukin-6 stimulation: contributions of Src- and receptor-tyrosine kinases, *J Biol Chem* 280, 31478–31488 (2005)
- Starzynski RR, Lipinski P, Drapier JC, Diet A, Smuda E, Bartlomieiczuk T, Gralak MA, and Kruszewski M, Down-regulation of iron regulatory protein 1 activities and expression in superoxide dismutase 1 knock-out mice is not associated with alterations in iron metabolism, *J Biol Chem* 280, 4207–4212 (2005)
- Su X and Dowhan W, Translational regulation of nuclear gene *COX4* expression by mitochondrial content of phosphatidylglycerol and cardiolipin in *Saccharomyces cerevisiae*, *Mol Cell Biol* 26, 743–753 (2006)
- Subramanian T, Vijayalingam S, and Chinnadurai G, Genetic identification of adenovirus type 5 genes that influence viral spread, *J Virol* 80, 2000–2012 (2006)
- Wang Y and Li G, ING3 promotes UV-induced apoptosis via fas/caspase-8 pathway in melanoma cells, *J Biol Chem*, M511309200, published electronically March 6 (2006)
- Youn S, Collisson EW, and Machamer CE, Contribution of trafficking signals in the cytoplasmic tail of the infectious bronchitis virus spike protein to virus infection, *J Virol* 79, 13209–13217 (2005)
- Zawadzka-Skomial J, Markiewicz Z, Nguyen-Disteche M, Devreese B, Frere JM, and Terrak M, Characterization of the bifunctional glycosyltransferase/acyltransferase penicillin-binding protein 4 of *Listeria monocytogenes*, *J Bacteriol* 188, 1875–1881 (2006)
- Zhou Q and Amar S, 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 (2006)