

The Discovery Series™ PDQuest™ 2-D Analysis Software Citations

Bulletin 2629 Rev C

- Acestor N, Masina S, Walker J, Saravia NG, Fasel N, and Quadroni M, Establishing two-dimensional gels for the analysis of *Leishmania* proteomes, *Proteomics* 2, 877–879 (2002)
- Ahmed N, Barker G, Oliva K, Garfin D, Talmadge K, Georgiou H, Quinn M, and Rice G, An approach to remove albumin for the proteomic analysis of low abundance biomarkers in human serum, *Proteomics* 3, 1980–1987 (2003)
- Ahmed N, Oliva K, Wang Y, Quinn M, and Rice G, Proteomic profiling of proteins associated with urokinase plasminogen activator receptor in a colon cancer cell line using an antisense approach, *Proteomics* 3, 288–298 (2003)
- Ahn WS, Kim KW, Bae SM, Yoon JH, Lee JM, Namkoong SE, Kim JH, Kim CK, Lee YJ, and Kim YW, Targeted cellular process profiling approach for uterine leiomyoma using cDNA microarray and gene ontology analysis, *Int J Exp Pathol* 84, 267–279 (2003)
- Alaiya A, Roblick U, Egevad L, Carlsson A, Franzen B, Volz D, Huwendiek S, Linder S, and Auer G, Polypeptide expression in prostate hyperplasia and prostate adenocarcinoma, *Anal Cell Pathol* 21, 1–9 (2000)
- Alaiya AA, Franzen B, Fujioka K, Moberger B, Schedvins K, Silfversvard C, Linder S, and Auer G, Phenotypic analysis of ovarian carcinoma: polypeptide expression in benign, borderline and malignant tumors, *Int J Cancer* 73, 678–683 (1997)
- Alaiya AA, Franzen B, Moberger B, Silfversward C, Linder S, and Auer G, Two-dimensional gel analysis of protein expression in ovarian tumors shows a low degree of intratumoral heterogeneity, *Electrophoresis* 20, 1039–1046 (1999)
- Alaiya AA, Oppermann M, Langridge J, Roblick U, Egevad L, Brindstedt S, Hellstrom M, Linder S, Bergman T, Jornvall H, and Auer G, Identification of proteins in human prostate tumor material by two-dimensional gel electrophoresis and mass spectrometry, *Cell Mol Life Sci* 58, 307–311 (2001)
- Alaiya AA, Roblick UJ, Franzen B, Bruch HP, and Auer G, Protein expression profiling in human lung, breast, bladder, renal, colorectal and ovarian cancers, *J Chromatogr B Analyt Technol Biomed Life Sci* 787, 207–222 (2003)
- Albers E, Laize V, Blomberg A, Hohmann S, and Gustafsson L, Ser3p (Yer081wp) and Ser33p (Yil074cp) are phosphoglycerate dehydrogenases in *Saccharomyces cerevisiae*, *J Biol Chem* 278, 10264–10272 (2003)
- Amrani M, Corbett J, Boateng SY, Dunn MJ, and Yacoub MH, Kinetics of induction and protective effect of heat-shock proteins after cardioplegic arrest, *Ann Thorac Surg* 61, 1407–1412 (1996)
- Antonucci F, Chilosi M, Parolini C, Hamdan M, Astner H, and Righetti PG, Two-dimensional molecular profiling of mantle cell lymphoma, *Electrophoresis* 24, 2376–2385 (2003)
- Axel H, Stephan MM, Thomas K, Peter A, Kai-Uwe K, and Sunita C, Multiple versus single ultrasound probes for tissue ablation by high intensity focused ultrasound: an experimental investigation, *J Urol* 169, 208 (2003)
- Aziz RK, Pabst MJ, Jeng A, Kansal R, Low DE, Nizet V, and Kotb M, Invasive M1T1 group A *Streptococcus* undergoes a phase-shift in vivo to prevent proteolytic degradation of multiple virulence factors by SpeB, *Mol Microbiol* 51, 123–134 (2004)
- Baik SC, Kim KM, Song SM, Kim DS, Jun JS, Lee SG, Song JY, Park JU, Kang HL, Lee WK, Cho MJ, Youn HS, Ko GH, and Rhee KH, Proteomic analysis of the sarcosine-inducible outer membrane fraction of *Helicobacter pylori* strain 26695, *J Bacteriol* 186, 949–955 (2004)
- Ballard DW, Bohnlein E, Lowenthal JW, Wano Y, Franza BR, and Greene WC, HTLV-I tax induces cellular proteins that activate the kappa B element in the IL-2 receptor alpha gene, *Science* 241, 1652–1655 (1988)
- Bardel J, Louwagie M, Jaquinod M, Jourdain A, Luche S, Rabilloud T, Macherel D, Garin J, and Bourguignon J, A survey of the plant mitochondrial proteome in relation to development, *Proteomics* 2, 880–898 (2002)
- Barker DL, Murray TF, Siebenaller JF, and Mpitsos GJ, Computer analyzed 2D gels reveal effects of muscarinic ligands on synthesis of specific proteins in the nervous system of *Pleurobranchaea californica*, *Soc Neurosci Abstr* 12, 400 (1986)
- Barnes S and Kim H, Nutriproteomics: identifying the molecular targets of nutritive and non-nutritive components of the diet, *J Biochem Mol Biol* 37, 59–74 (2004)
- Barry RC, Alsaker BL, Robison-Cox JF, and Dratz EA, Quantitative evaluation of sample application methods for semipreparative separations of basic proteins by two-dimensional gel electrophoresis, *Electrophoresis* 24, 3390–3404 (2003)

- Baty JW, Hampton MB, and Winterbourn CC, Detection of oxidant sensitive thiol proteins by fluorescence labeling and two-dimensional electrophoresis, *Proteomics* 2, 1261–1266 (2002)
- Bensmail H and Haoudi A, Postgenomics: proteomics and bioinformatics in cancer research, *J Biomed Biotechnol* 2003, 217–230 (2003)
- Berendt B, Detail and context in Web usage mining: coarsening and visualizing sequences, *WEBKDD 2001: Mining Web Log Data Across All Customer Touch Points* 2356, 1–24 (2002)
- Beresini MH and Epstein LB, Synergistic effects of interferon-gamma and tumor necrosis factor on the induction of specific polypeptides and the inhibition of growth of cervical carcinoma cells, *J Interferon Res* 7, 819 (1987)
- Beresini MH, Lempert MJ, and Epstein LB, Overlapping patterns of polypeptide induction by interleukin-1alpha (IL-1alpha), interleukin-1beta (IL-1beta), interferon-alpha (IFN-alpha), interferon-gamma (IFN-gamma), and tumor necrosis factor (TNF) in human fibroblasts, *Proc Amer Assoc Cancer Res* 28, 341 (1987)
- Beresini MH, Lempert MJ, and Epstein LB, Polypeptide induction indicates potential biochemical pathways in common in the mechanisms of action of IFN-alpha, IFN-gamma, IL-1 and TNF, *J Interferon Res* 7, 819 (1987)
- Beresini MH, Lempert MJ, and Epstein LB, Overlapping polypeptide induction in human fibroblasts in response to treatment with interferon-alpha, interferon-gamma, interleukin 1 alpha, interleukin 1 beta, and tumor necrosis factor, *J Immunol* 140, 485–493 (1988)
- Beresini MH, Sugarman BJ, Shepard HM, and Epstein LB, Synergistic induction of polypeptides by tumor necrosis factor and interferon-gamma in cells sensitive or resistant to tumor necrosis factor: assessment by computer based analysis of two-dimensional gels using the PDQUEST system, *Electrophoresis* 11, 232–241 (1990)
- Bershadsky AD, Gluck U, Denisenko ON, Sklyarova TV, Spector I, and Ben-Ze'ev A, The state of actin assembly regulates actin and vinculin expression by a feedback loop, *J Cell Sci* 108, 1183–1193 (1995)
- Binz PA, Hochstrasser DF, and Appel RD, Mass spectrometry-based proteomics: current status and potential use in clinical chemistry, *Clin Chem Lab Med* 41, 1540–1551 (2003)
- Blomberg A, Blomberg L, Norbeck J, Fey SJ, Larsen PM, Larsen M, Roepstorff P, Degand H, Boutry M, Posch A et al., Interlaboratory reproducibility of yeast protein patterns analyzed by immobilized pH gradient two-dimensional gel electrophoresis, *Electrophoresis* 16, 1935–1945 (1995)
- Blose SH, Cellular protein databases: a new approach to the study of genomic organization and function of the cell, *Leukemia* 2, 559 (1988)
- Blose SH, The use of computerized two-dimensional gels in characterizing hybridoma cell lines and monoclonal antibodies: applications in quality control and patent protection, *Hybridoma* 4, 66 (1985)
- Bohnlein E, Lowenthal JW, Siekevitz M, Ballard DW, Franza BR, and Greene WC, The same inducible nuclear proteins regulates mitogen activation of both the interleukin-2 receptor-alpha gene and type 1 HIV, *Cell* 53, 827–836 (1988)
- Bouchal P and Kucera I, Examination of membrane protein expression in *Paracoccus denitrificans* by two-dimensional gel electrophoresis, *J Basic Microbiol* 44, 17–22 (2004)
- Boxman IL, Hensbergen PJ, Van Der Schors RC, Bruynzeel DP, Tensen CP, and Ponc M, Proteomic analysis of skin irritation reveals the induction of HSP27 by sodium lauryl sulphate in human skin, *Br J Dermatol* 146, 777–785 (2002)
- Bradley BP, Shrader EA, Kimmel DG, and Meiller JC, Protein expression signatures: an application of proteomics, *Mar Environ Res* 54, 373–377 (2002)
- Broeckeaert F, Clippe A, Wattiez R, Falmagne P, and Bernard A, Lung hyperpermeability, Clara-cell secretory protein (CC16), and susceptibility to ozone of five inbred strains of mice, *Inhal Toxicol* 15, 1209–1230 (2003)
- Brophy B, Smolenski G, Wheeler T, Wells D, L'Huillier P, and Laible G, Cloned transgenic cattle produce milk with higher levels of beta-casein and kappa-casein, *Nat Biotechnol* 21, 157–162 (2003)
- Bruno ME, Borchers CH, Dial JM, Walker NJ, Hartis JE, Wetmore BA, Carl Barrett JC, Tomer KB, and Merrick BA, Effects of TCDD upon I kappa B and IKK subunits localized in microsomes by proteomics, *Arch Biochem Biophys* 406, 153–164 (2002)
- Caldini A, Moneti G, Fanelli A, Bruschettoni A, Mercurio S, Pieraccini G, Cini E, Ognibene A, Luceri F, and Messeri G, Epoetin alpha, epoetin beta and darbepoetin alfa: two-dimensional gel electrophoresis isoforms characterization and mass spectrometry analysis, *Proteomics* 3, 937–941 (2003)
- Calikowski TT, Meulia T, and Meier I, A proteomic study of the *Arabidopsis* nuclear matrix, *J Cell Biochem* 90, 361–378 (2003)
- Carboni L, Piubelli C, Righetti PG, Jansson B, and Domenici E, Proteomic analysis of rat brain tissue: Comparison of protocols for two-dimensional gel electrophoresis analysis based on different solubilizing agents, *Electrophoresis* 23, 4132–4141 (2002)
- Castellucci VF, Kandel ER, Kennedy TE, and Goelet P, Learning associated proteins (LAPs) following long-term sensitization of the gill and siphon withdrawal reflex in *Aplysia*, *Soc Neurosci Abstr* 13, 390 (1987)
- Castellucci VF, Kennedy TE, Kandel ER, and Goelet P, A quantitative analysis of 2-D gels identifies proteins in which labeling is increased following long-term sensitization in *Aplysia*, *Neuron* 1, 321–328 (1988)
- Cecconi D, Astner H, Donadelli M, Palmieri M, Missiaglia E, Hamdan M, Scarpa A, and Righetti PG, Proteomic analysis of pancreatic ductal carcinoma cells treated with 5-aza-2'-deoxycytidine, *Electrophoresis* 24, 4291–4303 (2003)

- Cecconi D, Scarpa A, Donadelli M, Palmieri M, Hamdan M, Astner H, and Righetti PG, Proteomic profiling of pancreatic ductal carcinoma cell lines treated with trichostatin-A, *Electrophoresis* 24, 1871–1878 (2003)
- Celis JE, Cruger D, Kiil J, Dejgaard K, Lauridsen JB, Ratz GP, Basse B, Celis A, Rasmussen HH, Bauw G et al., A two-dimensional gel protein database of noncultured total normal human epidermal keratinocytes: identification of proteins strongly up-regulated in psoriatic epidermis, *Electrophoresis* 11, 242–254 (1990)
- Celis JE, Cruger D, Kiil J, Lauridsen JB, Ratz G, Basse B, and Celis A, Identification of a group of proteins that are strongly up-regulated in total epidermal keratinocytes from psoriatic skin, *FEBS Lett* 262, 159–164 (1990)
- Celis JE, Dejgaard K, Madsen P, Leffers H, Gesser B, Honore B, Rasmussen HH, Olsen E, Lauridsen JB, Ratz G et al., The MRC-5 human embryonic lung fibroblast two-dimensional gel cellular protein database: quantitative identification of polypeptides whose relative abundance differs between quiescent, proliferating and SV40 transformed cells, *Electrophoresis* 11, 1072–1113 (1990)
- Celis JE, Gesser B, Dejgaard K, Honore B, Leffers H, Madsen P, Andersen A, Basse B, Celis A, Lauridsen JB et al., Two dimensional gel human protein databases offer a systematic approach to the study of cell proliferation and differentiation, *Int J Dev Biol* 33, 407–416 (1989)
- Celis JE, Gesser B, Rasmussen HH, Madsen P, Leffers H, Dejgaard K, Honore B, Olsen E, Ratz G, Lauridsen JB et al., Comprehensive two-dimensional gel protein databases offer a global approach to the analysis of human cells: the transformed amnion cells (AMA) master database and its link to genome DNA sequence data, *Electrophoresis* 11, 989–1071 (1990)
- Celis JE, Honore B, Bauw G and Vandekerckhove J, Comprehensive computerized 2D gel protein databases offer a global approach to the study of the mammalian cell, *Bioessays* 12, 93–97 (1990)
- Celis JE, Leffers H, Rasmussen HH, Madsen P, Honore B, Gesser B, Dejgaard K, Olsen E, Ratz GP, Lauridsen JB et al., The master two-dimensional gel database of human AMA cell proteins: towards linking protein and genome sequence and mapping information (update 1991), *Electrophoresis* 12, 765–801 (1991)
- Celis JE, Madsen P, Rasmussen HH, Leffers H, Honore B, Gesser B, Dejgaard K, Olsen E, Magnusson N, Kiil J et al., A comprehensive two-dimensional gel protein database of noncultured unfractionated normal human epidermal keratinocytes: towards an integrated approach to the study of cell proliferation, differentiation and skin diseases, *Electrophoresis* 12, 802–872 (1991)
- Celis JE, Ratz GP, and Celis A, Secreted proteins from normal and SV40 transformed human MRC-5 fibroblasts: toward establishing a database of human secreted proteins, *Leukemia* 1, 707–717 (1987)
- Celis JE, Ratz GP, Celis A, Madsen P, Gesser B, Kwee S, Madsen PS, Nielsen HV, Yde H, Lauridsen JB et al., Towards establishing comprehensive databases of cellular proteins from transformed human epithelial amnion cells (AMA) and normal peripheral blood mononuclear cells, *Leukemia* 2, 561–601 (1988)
- Celis JE, Ratz GP, Madsen P, Gesser B, Lauridsen JB, Hansen KP, Kwee S, Rasmussen HH, Nielsen HV, Cruger D et al., Computerized, comprehensive databases of cellular and secreted proteins from normal human embryonic lung MRC-5 fibroblasts: identification of transformation and/or proliferation sensitive proteins, *Electrophoresis* 10, 76–115 (1989)
- Celis JE, Ratz GP, Madsen P, Gesser B, Lauridsen JB, Kwee S, Rasmussen HH, Nielsen HV, Cruger D, Basse B et al., Comprehensive, human cellular protein databases and their implication for the study of genome organization and function, *FEBS Lett* 244, 247–254 (1989)
- Chew CS and Brown MR, Histamine increases phosphorylation of 27- and 40-kDa parietal cell proteins, *Am J Physiol* 253, G823–G829 (1987)
- Choe LH and Lee KH, Quantitative and qualitative measure of intralaboratory two-dimensional protein gel reproducibility and the effects of sample preparation, sample load, and image analysis, *Electrophoresis* 24, 3500–3507 (2003)
- Choi W, Yoo YJ, Kim M, Shin D, Jeon HB, and Choi W, Identification of proteins highly expressed in the hyphae of *Candida albicans* by two-dimensional electrophoresis, *Yeast* 20, 1053–1060 (2003)
- Chow KT, Pope MK, and Davies J, Characterization of a vanillic acid non-oxidative decarboxylation gene cluster from *Streptomyces* sp. D7, *Microbiology* 145, 2393–2403 (1999)
- Clack JW, Juhl M, Rice CA, Li J, and Witzmann FA, Proteomic analysis of transducin beta-subunit structural heterogeneity, *Electrophoresis* 24, 3493–3499 (2003)
- Collins H, Najafi H, Buettger C, Rombeau J, Settle RG, and Matschinsky FM, Identification of glucose response proteins in two biological models of beta-cell adaptation to chronic high glucose exposure, *J Biol Chem* 267, 1357–1366 (1992)
- Collins PJ, Juhl C, and Lognonne JL, Image analysis of 2D gels: considerations and insights, *Cell Mol Biol (Noisy-le-grand)* 40, 77–83 (1994)
- Coppée JY, Auger S, Turlin E, Sekowska A, Le Caer JP, Labas V, Vagner V, Danchin A, and Martin-Verstraete I, Sulfur-limitation-regulated proteins in *Bacillus subtilis*: a two-dimensional gel electrophoresis study, *Microbiology* 147, 1631–1640 (2001)
- Corbett JM, Dunn MJ, Posch A, and Gorg A, Positional reproducibility of protein spots in two-dimensional polyacrylamide gel electrophoresis using immobilised pH gradient isoelectric focusing in the first dimension: an interlaboratory comparison, *Electrophoresis* 15, 1205–1211 (1994)

- Corbett JM, Wheeler CH, and Dunn MJ, Coelectrophoresis of cardiac tissue from human, dog, rat and mouse: towards the establishment of an integrated two-dimensional protein database, *Electrophoresis* 16, 1524–1529 (1995)
- Cordwell SJ, Acquisition and archiving of information for bacterial proteomics: from sample preparation to database, *Methods Enzymol* 358, 207–227 (2002)
- Curran T, Rauscher FJ 3rd, Sambucetti LC, and Franza BR, Fos protein complex binds to the recognition sequence of the HeLa cell activator protein-1 (AP-1), in Franza BR, Cullen BR and Wong-Staal F (eds) *The Control of Human Retrovirus Gene Expression*, Cold Spring Harbor Press, New York (1988)
- Curran T, Sambucetti L, Rubino K, Franza BR, and Morgan JI, *c-Fos* induction: a possible nuclear switch in growth regulation, *ICSU Proceedings of the Miami Winter Symposium* (1986)
- Cutler P, Heald G, White IR, and Ruan J, A novel approach to spot detection for two-dimensional gel electrophoresis images using pixel value collection, *Proteomics* 3, 392–401 (2003)
- de Lamirande E and Gagnon C, Redox control of changes in protein sulfhydryl levels during human sperm capacitation, *Free Radic Biol Med* 35, 1271–1285 (2003)
- Decker ED, Zhang Y, Cocklin RR, Witzmann FA, and Wang M, Proteomic analysis of differential protein expression induced by ultraviolet light radiation in HeLa cells, *Proteomics* 3, 2019–2027 (2003)
- Desiderio DM and Zhan X, The human pituitary proteome: the characterization of differentially expressed proteins in an adenoma compared to a control, *Cell Mol Biol (Noisy-le-grand)* 49, 689–712 (2003)
- Diffie GM, Seversen EA, Stein TD, and Johnson JA, Microarray expression analysis of effects of exercise training: increase in atrial MLC-1 in rat ventricles, *Am J Physiol Heart Circ Physiol* 284, H830–H837 (2003)
- Ding SJ, Li Y, Tan YX, Jiang MR, Tian B, Liu YK, Shao XX, Ye SL, Wu JR, Zeng R, Wang HY, Tang ZY, and Xia QC, From proteomic analysis to clinical significance: overexpression of cytokeratin 19 correlates with hepatocellular carcinoma metastasis, *Mol Cell Proteomics* 3, 73–81 (2004)
- DiPaolo BR, Speicher KD, and Speicher DW, Identification of the amino acid mutations associated with human erythrocyte spectrin alpha II domain polymorphisms, *Blood* 82, 284–291 (1993)
- Dowsey AW, Dunn MJ, and Yang GZ, The role of bioinformatics in two-dimensional gel electrophoresis, *Proteomics* 3, 1567–1596 (2003)
- Drummel-Smith J, Brochu V, Girard I, Messier N, and Ouellette M, Proteome mapping of the protozoan parasite *Leishmania* and application to the study of drug targets and resistance mechanisms, *Mol Cell Proteomics* 2, 146–155 (2003)
- Duran MC, Mas S, Martin-Ventura JL, Meilhac O, Michel JB, Gallego-Delgado J, Lazaro A, Tunon J, Egido J, and Vivanco F, Proteomic analysis of human vessels: application to atherosclerotic plaques, *Proteomics* 3, 973–978 (2003)
- Dwek MV and Alaiya AA, Proteome analysis enables separate clustering of normal breast, benign breast and breast cancer tissues, *Br J Cancer* 89, 305–307 (2003)
- Dykes GA, Sampathkumar B, and Korber DR, Planktonic or biofilm growth affects survival, hydrophobicity and protein expression patterns of a pathogenic *Campylobacter jejuni* strain, *Int J Food Microbiol* 89, 1–10 (2003)
- Easty D, Hart IR, Patel K, Seymour C, Yacoub M, Domscheit A, Gunther S, Postel W, Gorg A, and Dunn MJ, Changes in protein expression during melanoma differentiation determined by computer analysis of 2-D gels, *Clin Exp Metastasis* 9, 221–230 (1991)
- Edvardsson U, Bergstrom M, Alexandersson M, Bamberg K, Ljung B, and Dahllöf B, Rosiglitazone (BRL49653), a PPARgamma-selective agonist, causes peroxisome proliferator-like liver effects in obese mice, *J Lipid Res* 40, 1177–1184 (1999)
- Edwards E, Kitt C, Oliver E, Finkelstein J, Wagster M, and McDonald WM, Depression and Parkinson's disease: a new look at an old problem, *Depress Anxiety* 16, 39–48 (2002)
- Eklund DM and Edqvist J, Localization of nonspecific lipid transfer proteins correlate with programmed cell death responses during endosperm degradation in *Euphorbia lagascae* seedlings, *Plant Physiol* 132, 1249–1259 (2003)
- El Fakhry Y, Ouellette M, and Papadopoulou B, A proteomic approach to identify developmentally regulated proteins in *Leishmania infantum*, *Proteomics* 2, 1007–1017 (2002)
- El Ghalbzouri A, Hensbergen P, Gibbs S, Kempenaar J, van der Schors R, and Ponc M, Fibroblasts facilitate re-epithelialization in wounded human skin equivalents, *Lab Invest* 84, 102–112 (2004)
- Encarnacion S, del Carmen Vargas M, Dunn MF, Davalos A, Mendoza G, Mora Y, and Mora J, *AniA* regulates reserve polymer accumulation and global protein expression in *Rhizobium etli*, *J Bacteriol* 184, 2287–2295 (2002)
- Encarnacion S, Guzman Y, Dunn MF, Hernandez M, del Carmen Vargas M, and Mora J, Proteome analysis of aerobic and fermentative metabolism in *Rhizobium etli* CE3, *Proteomics* 3, 1077–1085 (2003)
- Epstein LB, The odyssey of an interferonologist: from interferon-gamma to Down syndrome, tumor necrosis factor, and interleukin-1, and back again, *J Interferon Res* 7, 487–495 (1987)
- Eriksson P, Alipour H, Adler L, and Blomberg A, Rap1p-binding sites in the *Saccharomyces cerevisiae* *GPD1* promoter are involved in its response to NaCl, *J Biol Chem* 275, 29368–29376 (2000)
- Fazeli MS, Corbet J, Dunn MJ, Dolphin AC, and Bliss TV, Changes in protein synthesis accompanying long-term potentiation in the dentate gyrus in vivo, *J Neurosci* 13, 1346–1353 (1993)

- Fetterer RH and Barfield RC, Characterization of a developmentally regulated oocyst protein from *Eimeria tenella*, *J Parasitol* 89, 553–564 (2003)
- Florentz C, Molecular investigations on tRNAs involved in human mitochondrial disorders, *Biosci Rep* 22, 81–98 (2002)
- Franza BR Jr, Josephs SF, Gilman MZ, Ryan W, and Clarkson B, Characterization of cellular proteins recognizing the HIV enhancer using a microscale DNA-affinity precipitation assay, *Nature* 330, 391–395 (1987)
- Franza BR Jr, Maruyama K, Garrels JI, and Ruley HE, In vitro establishment is not a sufficient prerequisite for transformation by activated *ras* oncogenes, *Cell* 44, 409–418 (1986)
- Franza BR Jr, Rauscher FJ 3rd, Josephs SF, and Curran T, The Fos complex and Fos-related antigens recognize sequence elements that contain AP-1 binding sites, *Science* 239, 1150–1153 (1988)
- Franza BR Jr, Sambucetti LC, Cohen DR, and Curran T, Analysis of Fos protein complexes and Fos-related antigens by high-resolution two-dimensional gel electrophoresis, *Oncogene* 1, 213–221 (1987)
- Franzen B, Becker S, Mikkola R, Tidblad K, Tjernberg A, and Birnbaum S, Characterization of periplasmic *Escherichia coli* protein expression at high cell densities, *Electrophoresis* 20, 790–797 (1999)
- Franzen B, Duvefelt K, Jonsson C, Engelhardt B, Ottervald J, Wickman M, Yang Y, and Schuppe-Koistinen I, Gene and protein expression profiling of human cerebral endothelial cells activated with tumor necrosis factor- α , *Brain Res Mol Brain Res* 115, 130–146 (2003)
- Franzen B, Linder S, Alaiya AA, Eriksson E, Uruy K, Hirano T, Okuzawa K, and Auer G, Analysis of polypeptide expression in benign and malignant human breast lesions: down-regulation of cytokeratins, *Br J Cancer* 74, 1632–1638 (1996)
- Fujiwara M, Jin E, Ghazizadeh M, and Kawanami O, Differential expression of protease-activated receptors 1, 2, and 4 on human endothelial cells from different vascular sites, *Pathobiology* 71, 52–58 (2004)
- Gadgil HS, Pabst KM, Giorgianni F, Umstot ES, Desiderio DM, Beranova-Giorgianni S, Gerling IC, and Pabst MJ, Proteome of monocytes primed with lipopolysaccharide: analysis of the abundant proteins, *Proteomics* 3, 1767–1780 (2003)
- Galeva N and Altermann M, Comparison of one-dimensional and two-dimensional gel electrophoresis as a separation tool for proteomic analysis of rat liver microsomes: cytochromes P450 and other membrane proteins, *Proteomics* 2, 713–722 (2002)
- Gamble SC, Dunn MJ, Wheeler CH, Joiner MC, Adu-Poku A, and Arrand JE, Expression of proteins coincident with inducible radioprotection in human lung epithelial cells, *Cancer Res* 60, 2146–2151 (2000)
- Gao Z, Zuberi A, Quon MJ, Dong Z, and Ye J, Aspirin inhibits serine phosphorylation of insulin receptor substrate 1 in tumor necrosis factor-treated cells through targeting multiple serine kinases, *J Biol Chem* 278, 24944–24950 (2003)
- Garrels JI and Franza BR Jr, The REF52 protein database. Methods of database construction and analysis using the QUEST system and characterizations of protein patterns from proliferating and quiescent REF52 cells, *J Biol Chem* 264, 5283–5298 (1989)
- Garrels JI and Franza BR Jr, Transformation-sensitive and growth-related changes of protein synthesis in REF52 cells. A two-dimensional gel analysis of SV40-, adenovirus-, and Kirsten murine sarcoma virus-transformed rat cells using the REF52 protein database, *J Biol Chem* 264, 5299–5312 (1989)
- Garrels JI, Changes in protein synthesis during myogenesis in a clonal cell line, *Dev Biol* 73, 134–152 (1979)
- Garrels JI, Two dimensional gel electrophoresis and computer analysis of proteins synthesized by clonal cell lines, *J Biol Chem* 254, 7961–7977 (1979)
- Garrels JI, Quantitative two-dimensional gel electrophoresis of proteins, *Methods Enzymol* 100, 411–423 (1983)
- Garrels JI, The QUEST system for quantitative analysis of two-dimensional gels, *J Biol Chem* 264, 5269–5282 (1989)
- Garrels JI and Gibson W, Identification and characterization of multiple forms of actin, *Cell* 9, 793–805 (1976)
- Garrels JI and Shubert D, Modulation of protein synthesis by nerve growth factor, *J Biol Chem* 254, 7978–7985 (1979)
- Ghafouri B, Tagesson C, and Lindahl M, Mapping of proteins in human saliva using two-dimensional gel electrophoresis and peptide mass fingerprinting, *Proteomics* 3, 1003–1015 (2003)
- Gianazza E, Veber D, Eberini I, Buccellato FR, Mutti E, Sironi L, and Scalabrino G, Cobalamin (vitamin B12)-deficiency-induced changes in the proteome of rat cerebrospinal fluid, *Biochem J* 374, 239–246 (2003)
- Gmuender H, Kuratli K, Di Padova K, Gray CP, Keck W, and Evers S, Gene expression changes triggered by exposure of *Haemophilus influenzae* to novobiocin or ciprofloxacin: combined transcription and translation analysis, *Genome Res* 11, 28–42 (2001)
- Goldman MB, Knovich MA, and Goldman JN, Immunologic control of C3 gene expression in tissue macrophages, *J Immunol* 147, 4248–4255 (1991)
- Gonzalez-Barderas M, Gallego-Delgado J, Mas S, Duran MC, Lazaro A, Hernandez-Merida S, Egido J, and Vivanco F, Isolation of circulating human monocytes with high purity for proteomic analysis, *Proteomics* 4, 432–437 (2004)
- Gonzalez-Camacho F and Medina FJ, Identification of specific plant nucleolar phosphoproteins in a functional proteomic analysis, *Proteomics* 4, 407–417 (2004)
- Goodman R and Henderson AS, Exposure of salivary gland cells to low-frequency electromagnetic fields alters polypeptide synthesis, *Proc Natl Acad Sci USA* 85, 3928–3932 (1988)

- Govorun VM and Archakov AI, Proteomic technologies in modern biomedical science, *Biochemistry (Mosc)* 67, 1109–1123 (2002)
- Greene ND, Leung KY, Wait R, Begum S, Dunn MJ, and Copp AJ, Differential protein expression at the stage of neural tube closure in the mouse embryo, *J Biol Chem* 277, 41645–41651 (2002)
- Groblewski GE, Wishart MJ, Yoshida M, and Williams JA, Purification and identification of a 28-kDa calcium-regulated heat-stable protein. A novel secretagogue-regulated phosphoprotein in exocrine pancreas, *J Biol Chem* 271, 31502–31507 (1996)
- Gromov PS, Østergaard M, Gromova I, and Celis JE, Human proteomic databases: a powerful resource for functional genomics in health and disease, *Prog Biophys Mol Biol* 80, 3–22 (2002)
- Grünenfelder B, Rummel G, Vohradsky J, Röder D, Langen H, and Jenal U, Proteomic analysis of the bacterial cell cycle, *Proc Natl Acad Sci USA* 98, 4681–4686 (2001)
- Gu X, Meer SG, Miyagi M, Rayborn ME, Hollyfield JG, Crabb JW, and Salomon RG, Carboxyethylpyrrole protein adducts and autoantibodies, biomarkers for age-related macular degeneration, *J Biol Chem* 278, 42027–42035 (2003)
- Guedes Sde M, Vitorino R, Tomer K, Domingues MR, Correia AJ, Amado F, and Domingues P, *Drosophila melanogaster* larval hemolymph protein mapping, *Biochem Biophys Res Commun* 312, 545–554 (2003)
- Gustafsson JS, Blomberg A, and Rudemo M, Warping two-dimensional electrophoresis gel images to correct for geometric distortions of the spot pattern, *Electrophoresis* 23, 1731–1744 (2002)
- Ha GH, Lee SU, Kang DG, Ha NY, Kim SH, Kim J, Bae JM, Kim JW, and Lee CW, Proteome analysis of human stomach tissue: separation of soluble proteins by two-dimensional polyacrylamide gel electrophoresis and identification by mass spectrometry, *Electrophoresis* 23, 2513–2524 (2002)
- Hagell P, Whalley D, McKenna SP, and Lindvall O, Health status measurement in Parkinson's disease: validity of the PDQ-39 and Nottingham Health Profile, *Mov Disord* 18, 773–783 (2003)
- Hamdan M and Righetti PG, Assessment of protein expression by means of 2-D gel electrophoresis with and without mass spectrometry, *Mass Spectrom Rev* 22, 272–284 (2003)
- Harlow E, Franza BR Jr, and Schley C, Monoclonal antibodies specific for adenovirus early region 1A proteins: extensive heterogeneity in early region 1A products, *J Virol* 55, 533–546 (1985)
- Harlow E, Whyte P, Franza BR Jr, and Schley C, Association of adenovirus early-region 1A proteins with cellular polypeptides, *Mol Cell Biol* 6, 1579–1589 (1986)
- Harris CA, Derbin KS, Hunte-McDonough B, Krauss MR, Chen KT, Smith DM, and Epstein LB, Manganese superoxide dismutase is induced by IFN-gamma in multiple cell types. Synergistic induction by IFN-gamma and tumor necrosis factor or IL-1, *J Immunol* 147, 149–154 (1991)
- Hatada I, Hayashizaki Y, Hirotsune S, Komatsubara H, and Mukai T, A genomic scanning method for higher organisms using restriction sites as landmarks, *Proc Natl Acad Sci USA* 88, 9523–9527 (1991)
- Häusler KD, Quinn JMW, Horwood NJ, Ellis J, Lengel C, Martin TJ, Rubin JS, and Gillespie MT, Secreted frizzled-related protein (sFRP-1) inhibits TNFalpha- or RANKL-dependent osteoclast formation, *J Bone Miner Res* 17, S349 (2002)
- Hecht NB, Kleene KC, Distel RJ, and Silver LM, The differential expression of the actins and tubulins during spermatogenesis in the mouse, *Exp Cell Res* 153, 275–280 (1984)
- Hedegaard CJ, Thogersen IB, Enghild JJ, Klintworth GK, and Moller-Pedersen T, Transforming growth factor beta induced protein accumulation in granular corneal dystrophy type III (Reis-Bucklers dystrophy). Identification by mass spectrometry in 15 year old two-dimensional protein gels, *Mol Vis* 9, 355–359 (2003)
- Hermans C, Libotte V, Robin M, Clippe A, Wattiez R, Falmagne P, Langhendries JP, and Bernard A, Maternal tobacco smoking and lung epithelium-specific proteins in amniotic fluid, *Pediatr Res* 50, 487–494 (2001)
- Hernandez G, Lalioti V, Vandekerckhove J, Sierra JM, and Santaren JF, Identification and characterization of the expression of the translation initiation factor 4A (eIF4A) from *Drosophila melanogaster*, *Proteomics* 4, 316–326 (2004)
- Higginbotham JW, Smith JS, and Smith OS, Quantitative analysis of two-dimensional protein profiles of inbred lines of maize (*Zea mays* L.), *Electrophoresis* 12, 425–431 (1991)
- Hinerfeld D, Traini MD, Weinberger RP, Cochran B, Doctrow SR, Harry J, and Melov S, Endogenous mitochondrial oxidative stress: neurodegeneration, proteomic analysis, specific respiratory chain defects, and efficacious antioxidant therapy in superoxide dismutase 2 null mice, *J Neurochem* 88, 657–667 (2004)
- Hobbs JA, May R, Tanousis K, McNeill E, Mathies M, Gebhardt C, Henderson R, Robinson MJ, and Hogg N, Myeloid cell function in MRP-14 (S100A9) null mice, *Mol Cell Biol* 23, 2564–2576 (2003)
- Horiguchi A, Oya M, Uchida A, Marumo K, and Murai M, Elevated Akt activation and its impact on clinicopathological features of renal cell carcinoma, *J Urol* 169, 710–713 (2003)
- Houtman R, Krijgsveld J, Kool M, Romijn EP, Redegeld FA, Nijkamp FP, Heck AJ, and Humphery-Smith I, Lung proteome alterations in a mouse model for nonallergic asthma, *Proteomics* 3, 2008–2018 (2003)
- Hu L, Evers S, Lu ZH, Shen Y, and Chen J, Two-dimensional protein database of human pancreas, *Electrophoresis* 25, 512–518 (2004)

- Huang L, Li B, Luo C, Xie J, Chen P, and Liang S, Proteome comparative analysis of gynogenetic haploid and diploid embryos of goldfish (*Carassius auratus*), *Proteomics* 4, 235–243 (2004)
- Hunter T and Garrels JI, Characterization of the mRNAs for alpha-, beta- and gamma-actin, *Cell* 12, 767–781 (1977)
- Ishikawa N, Taniguchi-Seto H, Munakata Y, Takagi Y, Shimada N, and Kimura N, Multiple transcripts for rat nucleoside diphosphate kinase alpha isoform are structurally categorized into two groups that exhibit cell-specific expression and distinct translation potential, *J Biol Chem* 272, 3289–3295 (1997)
- Islam N, Tsujimoto H, and Hirano H, Wheat proteomics: relationship between fine chromosome deletion and protein expression, *Proteomics* 3, 307–316 (2003)
- Islam N, Tsujimoto H, and Hirano H, Proteome analysis of diploid, tetraploid and hexaploid wheat: Towards understanding genome interaction in protein expression, *Proteomics* 3, 549–557 (2003)
- Islam N, Woo SH, Tsujimoto H, Kawasaki H, and Hirano H, Proteome approaches to characterize seed storage proteins related to ditelocentric chromosomes in common wheat (*Triticum aestivum* L.), *Proteomics* 2, 1146–1155 (2002)
- Jackson HJ and Burgess PM, Personality disorders in the community: results from the Australian National Survey of Mental Health and Wellbeing Part II. Relationships between personality disorder, Axis I mental disorders and physical conditions with disability and health consultations, *Soc Psychiatry Psychiatr Epidemiol* 37, 251–260 (2002)
- Javor GT, Thiol-sensitive genes of *Escherichia coli*, *J Bacteriol* 171, 5607–5613 (1989)
- Jiang D, Ying W, Lu Y, Wan J, Zhai Y, Liu W, Zhu Y, Qiu Z, Qian X, and He F, Identification of metastasis-associated proteins by proteomic analysis and functional exploration of interleukin-18 in metastasis, *Proteomics* 3, 724–737 (2003)
- Jiang M, Jia L, Jiang W, Hu X, Zhou H, Gao X, Lu Z, and Zhang Z, Protein dysregulation in red blood cell membranes of type 2 diabetic patients, *Biochem Biophys Res Commun* 309, 196–200 (2003)
- Jin SH, Cho EH, Ko JE, Jung EH, Ahn B, Hahn JR, Kim JW, Kim CW, and Kim DR, Comparative analysis of nuclear proteins of B cells in different developmental stages, *Proteomics* 3, 2428–2436 (2003)
- Kennedy TE, Gawinowicz MA, Barzilai A, Kandel ER, and Sweatt JD, Sequencing of proteins from two-dimensional gels by using in situ digestion and transfer of peptides to polyvinylidene difluoride membranes: application to proteins associated with sensitization in *Aplysia*, *Proc Natl Acad Sci USA* 85, 7008–7012 (1988)
- Kennedy TE, Wager-Smith K, Barzilai A, Kandel ER, and Sweatt JD, Sequencing proteins from acrylamide gels, *Nature* 336, 499–500 (1988)
- Kernec F, Unlu M, Labeikovsky W, Minden JS, and Koretsky AP, Changes in the mitochondrial proteome from mouse hearts deficient in creatine kinase, *Physiol Genomics* 6, 117–128 (2001)
- Kim H, Page GP, and Barnes S, Proteomics and mass spectrometry in nutrition research, *Nutrition* 20, 155–165 (2004)
- Kim IY, Ahn HJ, Zelner DJ, Park L, Sensibar JA, and Lee C, Expression and localization of transforming growth factor-beta receptors type I and type II in the rat ventral prostate during regression, *Mol Endocrinol* 10, 107–115 (1996)
- Kim J, Kim SH, Lee SU, Ha GH, Kang DG, Ha NY, Ahn JS, Cho HY, Kang SJ, Lee YJ, Hong SC, Ha WS, Bae JM, Lee CW, and Kim JW, Proteome analysis of human liver tumor tissue by two-dimensional gel electrophoresis and matrix assisted laser desorption/ionization-mass spectrometry for identification of disease-related proteins, *Electrophoresis* 23, 4142–4156 (2002)
- Kim SK, Yoo JI, Cho BK, Hong SJ, Kim YK, Moon JA, Kim JH, Chung YN, and Wang KC, Elevation of CRABP-I in the cerebrospinal fluid of patients with Moyamoya disease, *Stroke* 34, 2835–2841 (2003)
- Kjaersgard IV and Jessen F, Proteome analysis elucidating post-mortem changes in cod (*Gadus morhua*) muscle proteins, *J Agric Food Chem* 51, 3985–3991 (2003)
- Knecht M, Regitz-Zagrosek V, Pleissner KP, Emig S, Jungblut P, Hildebrandt A, and Fleck E, Dilated cardiomyopathy: computer-assisted analysis of endomyocardial biopsy protein patterns by two-dimensional gel electrophoresis, *Eur J Clin Chem Clin Biochem* 32, 615–624 (1994)
- Knecht M, Regitz-Zagrosek V, Pleissner KP, Jungblut P, Steffen C, Hildebrandt A, and Fleck E, Characterization of myocardial protein composition in dilated cardiomyopathy by two-dimensional gel electrophoresis, *Eur Heart J* 15D, 37–44 (1994)
- Knecht M, Shinohara O, and Catt KJ, Biosynthesis of cellular and secreted proteins during follicle-stimulating hormone-induced granulosa cell differentiation, *Endocrinology* 119, 1388–1396 (1986)
- Konig S, Schmidt O, Rose K, Thanos S, Besselmann M, and Zeller M, Sodium dodecyl sulfate versus acid-labile surfactant gel electrophoresis: Comparative proteomic studies on rat retina and mouse brain, *Electrophoresis* 24, 751–756 (2003)
- Korke R, Gatti Mde L, Lau AL, Lim JW, Seow, TK, Chung MC, and Hu WS, Large scale gene expression profiling of metabolic shift of mammalian cells in culture, *J Biotechnol* 107, 1–17 (2004)
- Krah A, Miehlke S, Pleissner KP, Zimny-Arndt U, Kirsch C, Lehn N, Meyer TF, Jungblut PR, and Aebischer T, Identification of candidate antigens for serologic detection of *Helicobacter pylori*-infected patients with gastric carcinoma, *Int J Cancer* 108, 456–463 (2004)
- Krauss MR and Collins PJ, Protein database development: identification of glycosylated and phosphorylated proteins in unfractionated rat fibroblast lysates, *Electrophoresis* 10, 158–163 (1989)

- Krauss MR and Collins PJ, The use of powerful software for the quantitation of two-dimensional gel electrophoresis data, *Am Biotechnol Lab* 8, 34–38 (1990)
- Krauss MR, Collins PJ, and Bloise SH, Computer-analysed 2-D electrophoresis, *Nature* 337, 669–670 (1989)
- Krauss MR, Collins PJ, and Bloise SH, A two-dimensional acrylamide gel electrophoresis/computer software approach to decoding the human genome, *Biotechniques* 8, 218–223 (1990)
- Krin E, Sismeiro O, Danchin A, and Bertin PN, The regulation of Enzyme IIA(Glc) expression controls adenylate cyclase activity in *Escherichia coli*, *Microbiology* 148, 1553–1559 (2002)
- Kuerer HM, Goldknopf IL, Fritsche H, Krishnamurthy S, Sheta EA, and Hunt KK, Identification of distinct protein expression patterns in bilateral matched pair breast ductal fluid specimens from women with unilateral invasive breast carcinoma. High-throughput biomarker discovery, *Cancer* 95, 2276–2282 (2002)
- Kuncewicz T, Sheta EA, Goldknopf IL, and Kone B, Proteomic analysis of S-nitrosylated proteins in mesangial cells, *Mol Cell Proteomics* 2, 156–163 (2003)
- Lasbury ME, Durant PJ, and Lee CH, Life cycle stage-specific and encystment protein profiles in *Pneumocystis carinii*, *J Eukaryot Microbiol* 50, 636 (2003)
- Latham KE, Bautista FD, Hirao Y, O'Brien MJ, and Eppig JJ, Comparison of protein synthesis patterns in mouse cumulus cells and mural granulosa cells: effects of follicle-stimulating hormone and insulin on granulosa cell differentiation in vitro, *Biol Reprod* 61, 482–492 (1999)
- Lau AT, He QY, and Chiu JF, Proteomic technology and its biomedical applications, *Sheng Wu Hua Xue Yu Sheng Wu Wu Li Xue Bao (Shanghai)* 35, 965–975 (2003)
- Lee E, Jin Y, Hwang R, Rhee Y, and Lim S, Mass spectrometric proteome analysis of RANKL-induced protein expression in osteoclast differentiation, *J Bone Miner Res* 17, S349 (2002)
- Lee HJ, Park KJ, Lee AY, Park SG, Park BC, Lee KH, and Park SJ, Regulation of *fur* expression by RpoS and Fur in *Vibrio vulnificus*, *J Bacteriol* 185, 5891–5896 (2003)
- Lee MS, Kim KW, Kim YH, and Lee GM, Proteome analysis of antibody-expressing CHO cells in response to hyperosmotic pressure, *Biotechnol Prog* 19, 1734–1741 (2003)
- Lee SC, Yoon TG, Yoo YI, Bang YJ, Kim HY, Jeoung DI, and Kim HJ, Analysis of spinal cord proteome in the rats with mechanical allodynia after the spinal nerve injury, *Biotechnol Lett* 25, 2071–2078 (2003)
- Lefranc D, Almeras L, Dubucquoi S, de Seze J, Vermersch P, and Prin L, Distortion of the self-reactive IgG antibody repertoire in multiple sclerosis as a new diagnostic tool, *J Immunol* 172, 669–678 (2004)
- Leszczynski D, Nylund R, Joenvaara S, and Reivinen J, Applicability of discovery science approach to determine biological effects of mobile phone radiation, *Proteomics* 4, 426–431 (2004)
- Levy F, Bulet P, and Ehret-Sabatier L, Proteomic analysis of the systemic immune response of *Drosophila*, *Mol Cell Proteomics* 3, 156–166 (2004)
- Li C, Chen Z, Xiao Z, Wu X, Zhan X, Zhang X, Li M, Li J, Feng X, Liang S, Chen P, and Xie JY, Comparative proteomics analysis of human lung squamous carcinoma, *Biochem Biophys Res Commun* 309, 253–260 (2003)
- Li ZB, Lehar M, Braga N, Westra W, Liu LH, and Flint PW, Study of human laryngeal muscle protein using two-dimensional electrophoresis and mass spectrometry, *Proteomics* 3, 1325–1334 (2003)
- Liu B, Wang J, Yu H, Chung P, and Bringham F, Estrogen inhibits PTH-stimulated osteoclastogenesis through prevention of PTH-induced osteoprotegerin downregulation in mouse marrow stromal cells, *J Bone Miner Res* 17, S348–S349 (2002)
- Liu JW, Shen JJ, Tanzillo-Swarts A, Bhatia B, Maldonado CM, Person MD, Lau SS, and Tang DG, Annexin II expression is reduced or lost in prostate cancer cells and its re-expression inhibits prostate cancer cell migration, *Oncogene* 22, 1475–1485 (2003)
- Liu X, Wu Y, Zehner ZE, Jackson-Cook C, and Ware JL, Proteomic analysis of the tumorigenic human prostate cell line M12 after microcell-mediated transfer of chromosome 19 demonstrates reduction of vimentin, *Electrophoresis* 24, 3445–3453 (2003)
- Ljung B, Bamberg K, Dahllof B, Kjellstedt A, Oakes ND, Ostling J, Svensson L, and Camejo G, AZ 242, a novel PPARalpha/gamma agonist with beneficial effects on insulin resistance and carbohydrate and lipid metabolism in ob/ob mice and obese Zucker rats, *J Lipid Res* 43, 1855–1863 (2002)
- Lonosky PM, Zhang X, Honavar VG, Dobbs DL, Fu A, and Rodermerl SR, A proteomic analysis of maize chloroplast biogenesis, *Plant Physiol* 134, 560–574 (2004)
- Lu HS, Jones MD, Shieh JH, Mendiaz EA, Feng D, Watler P, Narhi LO, and Langley KE, Isolation and characterization of a disulfide-linked human stem cell factor dimer. Biochemical, biophysical, and biological comparison to the noncovalently held dimer, *J Biol Chem* 271, 11309–11316 (1996)
- Luria A, Vegelyte-Avery V, Stith B, Tsvetkova NM, Wolkers WF, Crowe JH, Tablin F, and Nuccitelli R, Detergent-free domain isolated from *Xenopus* egg plasma membrane with properties similar to those of detergent-resistant membranes, *Biochemistry* 41, 13189–13197 (2002)
- MacFarlane AJ, Burghardt KM, Kelly J, Simell T, Simell O, Altosaar I, and Scott FW, A type 1 diabetes-related protein from wheat (*Triticum aestivum*). cDNA clone of a wheat storage globulin, Glb1, linked to islet damage, *J Biol Chem* 278, 54–63 (2003)

- Madi A, Pusztahelyi T, Punyiczki M, and Fesus L, The biology of the post-genomic era: the proteomics, *Acta Biol Hung* 54, 1–14 (2003)
- Madsen PS, Hokland M, Ellegaard J, Hokland P, Ratz GP, Celis A, and Celis JE, Major proteins in normal human lymphocyte subpopulations separated by fluorescence-activated cell sorting and analyzed by two-dimensional gel electrophoresis, *Leukemia* 2, 602–615 (1988)
- Magnusson LU, Nystrom T, and Farewell A, Underproduction of sigma 70 mimics a stringent response. A proteome approach, *J Biol Chem* 278, 968–973 (2003)
- Malmstrom J, Tufvesson E, Lofdahl CG, Hansson L, Marko-Varga G, and Westergren-Thorsson G, Activation of platelet-derived growth factor pathway in human asthmatic pulmonary-derived mesenchymal cells, *Electrophoresis* 24, 276–285 (2003)
- Malmstrom L, Malmstrom J, Marko-Varga G, and Westergren-Thorsson G, Proteomic 2DE database for spot selection, automated annotation, and data analysis, *J Proteome Res* 1, 135–138 (2002)
- Marengo E, Robotti E, Righetti PG, and Antonucci F, New approach based on fuzzy logic and principal component analysis for the classification of two-dimensional maps in health and disease. Application to lymphomas, *J Chromatogr A* 1004, 13–28 (2003)
- Marko-Varga G, Berglund M, Malmstrom J, Lindberg H, and Fehniger TE, Targeting hepatocytes from liver tissue by laser capture microdissection and proteomics expression profiling, *Electrophoresis* 24, 3800–3805 (2003)
- Martinet W, Schrijvers DM, De Meyer GR, Herman AG, and Kockx MM, Western array analysis of human atherosclerotic plaques: down regulation of apoptosis-linked gene 2, *Cardiovasc Res* 60, 259–267 (2003)
- Martinez-Rodriguez R, Toledano A, Alvarez MI, Turegano L, Colman O, Roses P, Gomez de Segura I, and De Miguel E, Chronic nicotine administration increases NGF-like immunoreactivity in frontoparietal cerebral cortex, *J Neurosci Res* 73, 708–716 (2003)
- Mathews MB, Bernstein RM, Franza BR Jr, and Garrels JI, Identity of the proliferating cell nuclear antigen and cyclin, *Nature* 309, 374–376 (1984)
- Matousek P, Hodny Z, Svandova I, and Svoboda P, Different methods of membrane domains isolation result in similar 2-D distribution patterns of membrane domain proteins, *Biochem Cell Biol* 81, 365–372 (2003)
- McClelland CM and Gullick WJ, Identification of surrogate markers for determining drug activity using proteomics, *Biochem Soc Trans* 31, 1488–1490 (2003)
- McGregor E, Kempster L, Wait R, Gosling M, Dunn MJ, and Powell JT, F-actin capping (CapZ) and other contractile saphenous vein smooth muscle proteins are altered by hemodynamic stress: a proteomic approach, *Mol Cell Proteomics* 3, 115–124 (2004)
- McLaughlin PJ, Stack BC Jr, Levin RJ, Fedok F, and Zagon IS, Defects in the opioid growth factor receptor in human squamous cell carcinoma of the head and neck, *Cancer* 97, 1701–1710 (2003)
- Melin P, Schnurer J, and Wagner EG, Proteome analysis of *Aspergillus nidulans* reveals proteins associated with the response to the antibiotic concanamycin A, produced by *Streptomyces* species, *Mol Genet Genomics* 267, 695–702 (2002)
- Michishita E, Kurahashi T, Suzuki T, Fukuda M, Fujii M, Hirano H, and Ayusawa D, Changes in nuclear matrix proteins during the senescence-like phenomenon induced by 5-chloro-deoxyuridine in HeLa cells, *Exp Gerontol* 37, 885–890 (2002)
- Mikulik K, Bobek J, Bezouskova S, Benada O, and Kofronova O, Expression of proteins and protein kinase activity during germination of aerial spores of *Streptomyces granaticolor*, *Biochem Biophys Res Commun* 299, 335–342 (2002)
- Molin M, Larsson T, Karlsson KA, and Blomberg A, Fragmentation of dihydroxyacetone kinase 1 from *Saccharomyces cerevisiae* indicates a two-domain structure, *Proteomics* 3, 752–763 (2003)
- Moritz B and Meyer HE, Approaches for the quantification of protein concentration ratios, *Proteomics* 3, 2208–2220 (2003)
- Murray EJ, Murray SS, and Manolagas SC, Two-dimensional gel autoradiographic analyses of the effects of 1,25-dihydroxy-cholecalciferol on protein synthesis in clonal rat osteosarcoma cells, *Endocrinology* 126, 2679–2692 (1990)
- Neidhardt FC, Report of workshop on cellular protein databases derived from two-dimensional polyacrylamide gel electrophoresis, *Electrophoresis* 10, 73–75 (1989)
- Neidhardt FC, Appleby DB, Sankar P, Hutton ME, and Phillips TA, Genomically linked cellular protein databases derived from two-dimensional polyacrylamide gel electrophoresis, *Electrophoresis* 10, 116–122 (1989)
- Nicodeme M, Perrin C, Hols P, Bracquart P, and Gaillard JL, Identification of an iron-binding protein of the Dps family expressed by *Streptococcus thermophilus*, *Curr Microbiol* 48, 51–56 (2004)
- Nilsson A, Norbeck J, Oelz R, Blomberg A, and Gustafsson L, Fermentative capacity after cold storage of baker's yeast is dependent on the initial physiological state but not correlated to the levels of glycolytic enzymes, *Int J Food Microbiol* 71, 111–124 (2001)
- Nishigaki R, Shinohara T, Toda T, Omori A, Ichinose S, Itoh M, Shirayoshi Y, Kurimasa A, and Oshimura M, An extra human chromosome 21 reduces *mhc-2a* expression in chimeric mice and Down syndrome, *Biochem Biophys Res Commun* 295, 112–118 (2002)
- Nishihara JC and Champion KM, Quantitative evaluation of proteins in one- and two-dimensional polyacrylamide gels using a fluorescent stain, *Electrophoresis* 23, 2203–2215 (2002)

- Nishimura K, Arai Y, Nakai Y, Shimizu K, Masashi N, Inoue H, Tokizane T, Takaha N, Nomomura N, and Okuyama A, Effects of dexamethasone on renal cell carcinoma, *J Urol* 169, 208 (2003)
- Noel-Georis I, Vallaeyts T, Chauvaux R, Monchy S, Falmagne P, Mergeay M, and Wattiez R, Global analysis of the *Ralstonia metallidurans* proteome: prelude for the large-scale study of heavy metal response, *Proteomics* 4, 151–179 (2004)
- Norbeck J and Blomberg A, Metabolic and regulatory changes associated with growth of *Saccharomyces cerevisiae* in 1.4 M NaCl. Evidence for osmotic induction of glycerol dissimilation via the dihydroxyacetone pathway, *J Biol Chem* 272, 5544–5554 (1997)
- Norbeck J, Pahlman AK, Akhtar N, Blomberg A, and Adler L, Purification and characterization of two isoenzymes of DL-glycerol-3-phosphatase from *Saccharomyces cerevisiae*. Identification of the corresponding *GPP1* and *GPP2* genes and evidence for osmotic regulation of Gpp2p expression by the osmosensing mitogen-activated protein kinase signal transduction pathway, *J Biol Chem* 271, 13875–13881 (1996)
- Nouwens AS, Beatson SA, Whitchurch CB, Walsh BJ, Schweizer HP, Mattick JS, and Cordwell SJ, Proteome analysis of extracellular proteins regulated by the *las* and *rhl* quorum sensing systems in *Pseudomonas aeruginosa* PAO1, *Microbiology* 149, 1311–1322 (2003)
- Nouwens AS, Willcox MD, Walsh BJ, and Cordwell SJ, Proteomic comparison of membrane and extracellular proteins from invasive (PAO1) and cytotoxic (6206) strains of *Pseudomonas aeruginosa*, *Proteomics* 2, 1325–1346 (2002)
- Novotna J, Vohradsky J, Berndt P, Gramajo H, Langen H, Li XM, Minas W, Orsaria L, Roeder D, and Thompson CJ, Proteomic studies of diauxic lag in the differentiating prokaryote *Streptomyces coelicolor* reveal a regulatory network of stress-induced proteins and central metabolic enzymes, *Mol Microbiol* 48, 1289–1303 (2003)
- Obermeyer N, Janson N, Bergmann J, Buck F, and Ito WD, Proteome analysis of migrating versus nonmigrating rat heart endothelial cells reveals distinct expression patterns, *Endothelium* 10, 167–178 (2003)
- Oguri T, Takahata I, Katsuta K, Nomura E, Hidaka M, and Inagaki N, Proteome analysis of rat hippocampal neurons by multiple large gel two-dimensional electrophoresis, *Proteomics* 2, 666–672 (2002)
- Oppermann M, Cols N, Nyman T, Helin J, Saarinen J, Byman I, Toran N, Alaiya AA, Bergman T, Kalkkinen N, Gonzalez-Duarte R, and Jornvall H, Identification of foetal brain proteins by two-dimensional gel electrophoresis and mass spectrometry comparison of samples from individuals with or without chromosome 21 trisomy, *Eur J Biochem* 267, 4713–4719 (2000)
- Osaka H, Wang YL, Takada K, Takizawa S, Setsuie R, Li H, Sato Y, Nishikawa K, Sun YJ, Sakurai M, Harada T, Hara Y, Kimura I, Chiba S, Namikawa K, Kiyama H, Noda M, Aoki S, and Wada K, Ubiquitin carboxy-terminal hydrolase L1 binds to and stabilizes monoubiquitin in neuron, *Hum Mol Genet* 12, 1945–1958 (2003)
- Otto K, Norbeck J, Larsson T, Karlsson KA, and Hermansson M, Adhesion of type 1-fimbriated *Escherichia coli* to abiotic surfaces leads to altered composition of outer membrane proteins, *J Bacteriol* 183, 2445–2453 (2001)
- Ozdemir A, Bozkurt S, Akdas A, and Simsek F, Prognostic value and histopathological relation of micro vessel density, Vascular Endothelial Growth Factor and Thrombospondin-1 in renal cell carcinoma, *J Urol* 169, 208 (2003)
- Özkan M, Dilek FB, Yetis Ü, and Özcengiz G, Nutritional and cultural parameters influencing antidipteran delta-endotoxin production, *Res Microbiol* 154, 49–53 (2003)
- Pang JX, Ginanni N, Dongre AR, Hefta SA, and Opitek GJ, Biomarker discovery in urine by proteomics, *J Proteome Res* 1, 161–169 (2002)
- Park M, Kang H, Lim S, Lee CT, and Kim O, Differential protein expressions induced by adenovirus-mediated p16 gene transfer in Balb/c nude mouse, *Proteomics* 3, 2412–2419 (2003)
- Paulson L, Martin P, Persson A, Nilsson CL, Ljung E, Westman-Brinkmalm A, Eriksson PS, Blennow K, and Davidsson P, Comparative genome- and proteome analysis of cerebral cortex from MK-801-treated rats, *J Neurosci Res* 71, 526–533 (2003)
- Peng C, Liang SP, Tan C, Lv HB, Huang H, Zhou M, Wang R, Li XL, and Li GY, [Researching a novel NPC-related candidate suppressor gene BRD7 by two-dimensional gel electrophoresis and MALDI-TOF-MS], *Sheng Wu Hua Xue Yu Sheng Wu Wu Li Xue Bao (Shanghai)* 35, 816–822 (2003)
- Periago PM, Abee T, and Wouters JA, Analysis of the heat-adaptive response of psychrotrophic *Bacillus weihenstephanensis*, *Int J Food Microbiol* 79, 17–26 (2002)
- Periago PM, van Schaik W, Abee T, and Wouters JA, Identification of proteins involved in the heat stress response of *Bacillus cereus* ATCC 14579, *Appl Environ Microbiol* 68, 3486–3495 (2002)
- Pferdeort VA, Wood TK, and Reardon KF, Proteomic changes in *Escherichia coli* TG1 after metabolic engineering for enhanced trichloroethene biodegradation, *Proteomics* 3, 1066–1069 (2003)
- Phan-Thanh L, Proteomic analysis of response to acid in *Listeria monocytogenes*, *Methods Enzymol* 358, 256–276 (2002)
- Pitarch A, Sanchez M, Nombela C, and Gil C, Analysis of the *Candida albicans* proteome. I. Strategies and applications, *J Chromatogr B Analyt Technol Biomed Life Sci* 787, 101–128 (2003)

- Pitarch A, Sanchez M, Nombela C, and Gil C, Analysis of the *Candida albicans* proteome. II. Protein information technology on the Net (update 2002), *J Chromatogr B Analyt Technol Biomed Life Sci* 787, 129–148 (2003)
- Pleissner KP, Regitz-Zagrosek V, Weise C, Neuss M, Krudewagen B, Soding P, Buchner K, Hucho F, Hildebrandt A, and Fleck E, Chamber-specific expression of human myocardial proteins detected by two-dimensional gel electrophoresis, *Electrophoresis* 16, 841–850 (1995)
- Plymoth A, Lofdahl CG, Ekberg-Jansson A, Dahlback M, Lindberg H, Fehniger TE, and Marko-Varga G, Human bronchoalveolar lavage: Biofluid analysis with special emphasis on sample preparation, *Proteomics* 3, 962–972 (2003)
- Poland J, Bohme A, Schubert K, and Sinha P, Revisiting electroblotting of immobilized pH gradient gels: A new protocol for studying post-translational modification of proteins, *Electrophoresis* 23, 4067–4071 (2002)
- Poland J, Cahill MA, and Sinha P, Isoelectric focusing in long immobilized pH gradient gels to improve protein separation in proteomic analysis, *Electrophoresis* 24, 1271–1275 (2003)
- Poland J, Urbani A, Lage H, Schnolzer M, and Sinha P, Study of the development of thermoresistance in human pancreatic carcinoma cell lines using proteome analysis, *Electrophoresis* 25, 173–183 (2004)
- Polevoda B, Norbeck J, Takakura H, Blomberg A, and Sherman F, Identification and specificities of N-terminal acetyltransferases from *Saccharomyces cerevisiae*, *EMBO J* 18, 6155–6168 (1999)
- Pyo J, Hwang SI, Oh J, Lee SJ, Kang SC, Kim JS, and Lim J, Characterization of a bovine pregnancy-associated protein using two dimensional gel electrophoresis, N-terminal sequencing and mass spectrometry, *Proteomics* 3, 2420–2427 (2003)
- Rakwal R and Agrawal GK, Rice proteomics: Current status and future perspectives, *Electrophoresis* 24, 3378–3389 (2003)
- Ramachandran A, Ceasar E and Darley-Usmar VM, Chronic exposure to nitric oxide alters the free iron pool in endothelial cells: role of mitochondrial respiratory complexes and heat shock proteins, *Proc Natl Acad Sci USA* 101, 384–389 (2004)
- Rasmussen HH, Orntoft TF, Wolf H, and Celis JE, Towards a comprehensive database of proteins from the urine of patients with bladder cancer, *J Urol* 155, 2113–2119 (1996)
- Rauscher FJ 3rd, Cohen DR, Curran T, Bos TJ, Vogt PK, Bohmann D, Tjian R, and Franza BR Jr, Fos-associated protein p39 is the product of the *jun* proto-oncogene, *Science* 240, 1010–1016 (1988)
- Rautajoki K, Nyman TA, and Lahesmaa R, Proteome characterization of human T helper 1 and 2 cells, *Proteomics* 4, 84–92 (2004)
- Ray P, Chen L, Tyurin VA, Kagan VE, and Witzmann FA, Proteomic and inducible transgenic approaches to study disease processes, *Am J Respir Cell Mol Biol* 29, S42–S46 (2003)
- Ren Y and Smith A, Mechanism of metallothionein gene regulation by heme-hemopexin. Roles of protein kinase C, reactive oxygen species, and cis-acting elements, *J Biol Chem* 270, 23988–23995 (1995)
- Ribom D, Westman-Brinkmalm A, Smits A, and Davidsson P, Elevated levels of alpha-2-Heremans-Schmid glycoprotein in CSF of patients with low-grade gliomas, *Tumour Biol* 24, 94–99 (2003)
- Riffeser M and Hock B, Vitellogenin levels in mussel hemolymph — a suitable biomarker for the exposure to estrogens? *Comp Biochem Physiol C Toxicol Pharmacol* 132, 75–84 (2002)
- Righetti PG, Castagna A, Antonucci F, Piubelli C, Cecconi D, Campostrini N, Zanusso G, and Monaco S, The proteome: anno Domini 2002, *Clin Chem Lab Med* 41, 425–438 (2003)
- Rodriguez J, Agudo M, Van Damme J, Vandekerckhove J, and Santaren JF, Polypeptides differentially expressed in imaginal discs define the peroxiredoxin family of genes in *Drosophila*, *Eur J Biochem* 267, 487–497 (2000)
- Rogers M, Graham J and Tonge RP, Using statistical image models for objective evaluation of spot detection in two-dimensional gels, *Proteomics* 3, 879–886 (2003)
- Rosengren AT, Salmi JM, Aittokallio T, Westerholm J, Lahesmaa R, Nyman TA, and Nevalainen OS, Comparison of PDQuest and Progenesis software packages in the analysis of two-dimensional electrophoresis gels, *Proteomics* 3, 1936–1946 (2003)
- Rujoi M, Jin J, Borchman D, Tang D, and Yappert MC, Isolation and lipid characterization of cholesterol-enriched fractions in cortical and nuclear human lens fibers, *Invest Ophthalmol Vis Sci* 44, 1634–1642 (2003)
- Saintier D, Burde MA, Rey JM, Maudelonde T, de Vernejoul MC, and Cohen-Solal ME, Estrogen receptor beta expression increases with osteoclast differentiation but is not regulated by estradiol and RANK-L, *J Bone Miner Res* 17, S349 (2002)
- Saintier D, de Vernejoul MC, and Cohen-Solal ME, Estradiol down regulates beta3 integrin expression on differentiating and mature human osteoclasts, *J Bone Miner Res* 17, S349 (2002)
- Sakaguchi H, Miyagi M, Darrow RM, Crabb JS, Hollyfield JG, Organisciak DT, and Crabb JW, Intense light exposure changes the crystallin content in retina, *Exp Eye Res* 76, 131–133 (2003)
- Sakharov DV, Bunschoten A, van Weelden H, and Wirtz KW, Photodynamic treatment and H₂O₂-induced oxidative stress result indifferent patterns of cellular protein oxidation, *Eur J Biochem* 270, 4859–4865 (2003)
- Salmi J, Aittokallio T, Westerholm J, Griese M, Rosengren A, Nyman TA, Lahesmaa R, and Nevalainen O, Hierarchical grid transformation for image warping in the analysis of two-dimensional electrophoresis gels, *Proteomics* 2, 1504–1515 (2002)

- Salusjarvi L, Poutanen M, Pitkanen JP, Koivistoinen H, Aristidou A, Kalkkinen N, Ruohonen L, and Penttila M, Proteome analysis of recombinant xylose-fermenting *Saccharomyces cerevisiae*, *Yeast* 20, 295–314 (2003)
- Santamaria E, Avila MA, Latasa MU, Rubio A, Martin-Duce A, Lu SC, Mato JM, and Corrales FJ, Functional proteomics of nonalcoholic steatohepatitis: mitochondrial proteins as targets of S-adenosylmethionine, *Proc Natl Acad Sci USA* 100, 3065–3070 (2003)
- Sarry JE, Sommerer N, Sauvage FX, Bergoin A, Rossignol M, Albangac G, and Romieu C, Grape berry biochemistry revisited upon proteomic analysis of the mesocarp, *Proteomics* 4, 201–215 (2004)
- Sawicki G, Dakour J, and Morrish DW, Functional proteomics of neurokinin B in the placenta indicates a novel role in regulating cytotrophoblast antioxidant defences, *Proteomics* 3, 2044–2051 (2003)
- Schlags W, Lachmann B, Walther M, Kratzel M, and Noe CR, Two-dimensional electrophoresis of recombinant human erythropoietin: a future method for the European Pharmacopoeia? *Proteomics* 2, 679–682 (2002)
- Seike M, Kondo T, Mori Y, Gemma A, Kudoh S, Sakamoto M, Yamada T, and Hirohashi S, Proteomic analysis of intestinal epithelial cells expressing stabilized beta-catenin, *Cancer Res* 63, 4641–4647 (2003)
- Seliger B, Menig M, Lichtenfels R, Atkins D, Bukur J, Halder TM, Kersten M, Harder A, Ackermann A, Beck J, Muehlenweg B, Brenner W, Melchior S, Kellner R, and Lottspeich F, Identification of markers for the selection of patients undergoing renal cell carcinoma-specific immunotherapy, *Proteomics* 3, 979–990 (2003)
- Shefcheck K, Yao X, and Fenselau C, Fractionation of cytosolic proteins on an immobilized heparin column, *Anal Chem* 75, 1691–1698 (2003)
- Shekouh AR, Thompson CC, Prime W, Campbell F, Hamlett J, Herrington CS, Lemoine NR, Crnogorac-Jurcevic T, Buechler MW, Friess H, Neoptolemos JP, Pennington SR, and Costello E, Application of laser capture microdissection combined with two-dimensional electrophoresis for the discovery of differentially regulated proteins in pancreatic ductal adenocarcinoma, *Proteomics* 3, 1988–2001 (2003)
- Shen WP, Aldrich TH, Venta-Perez G, Franza BR Jr, and Furth ME, Expression of normal and mutant ras proteins in human acute leukemia, *Oncogene* 1, 157–165 (1987)
- Sheng H, Enghild JJ, Bowler R, Patel M, Batinic-Haberle I, Calvi CL, Day BJ, Pearlstein RD, Crapo JD, and Warner DS, Effects of metalloporphyrin catalytic antioxidants in experimental brain ischemia, *Free Radic Biol Med* 33, 947–961 (2002)
- Shi CZ, Collins HW, Buettger CW, Garside WT, Matschinsky FM, and Heyner S, Insulin family growth factors have specific effects on protein synthesis in preimplantation mouse embryos, *Mol Reprod Dev* 37, 398–406 (1994)
- Shi CZ, Collins HW, Garside WT, Buettger CW, Matschinsky FM, and Heyner S, Protein databases for compacted eight-cell and blastocyst-stage mouse embryos, *Mol Reprod Dev* 37, 34–47 (1994)
- Shi T, Liou LS, Dong F, and Didonato JA, Comparison of protein profiles in renal tissues and their primary cell lines, *J Urol* 169, 201 (2003)
- Shi T, Liou LS, Dong F, and Didonato JA, Protein alterations in human renal cell carcinoma analyzed by two-dimensional polyacrylamide gel electrophoresis with correlation to RNA expression, *J Urol* 169, 208 (2003)
- Shimazui T, Yoshikawa K, Uemura H, Hirao Y, Saga S, and Akaza H, Messenger RNA level of cadherin-6 in peripheral blood is associated with site of metastasis and occurrence of future metastasis in renal cell carcinoma, *J Urol* 169, 207–208 (2003)
- Silver LM, Uman J, Danska J, and Garrels JI, A diversified set of testicular cell proteins specified by genes within the mouse t complex, *Cell* 35, 35–45 (1983)
- Simpson DM, Beynon RJ, Robertson DH, Loughran MJ, and Haywood S, Copper-associated liver disease: a proteomics study of copper challenge in a sheep model, *Proteomics* 4, 524–536 (2004)
- Singh SK, Chhiber SS, Vohra H, and Sharma SK, Serum transforming growth factor beta 1 in patients of renal cell carcinoma, *J Urol* 169, 201 (2003)
- Sinha P, Hutter G, Kottgen E, Dietel M, Schadendorf D, and Lage H, Increased expression of annexin I and thioredoxin detected by two-dimensional gel electrophoresis of drug resistant human stomach cancer cells, *J Biochem Biophys Methods* 37, 105–116 (1998)
- Sinha P, Poland J, Kohl S, Schnolzer M, Helmbach H, Hutter G, Lage H, and Schadendorf D, Study of the development of chemoresistance in melanoma cell lines using proteome analysis, *Electrophoresis* 24, 2386–2404 (2003)
- Sironi L, Tremoli E, Miller I, Guerrini U, Calvio AM, Eberini I, Gemeiner M, Asdente M, Paoletti R, and Gianazza E, Acute-phase proteins before cerebral ischemia in stroke-prone rats: identification by proteomics, *Stroke* 32, 753–760 (2001)
- Smit E, Jager D, Martinez B, Tielen FJ, and Pouwels PH, Structural and functional analysis of the S-layer protein crystallisation domain of *Lactobacillus acidophilus* ATCC 4356: evidence for protein-protein interaction of two subdomains, *J Mol Biol* 324, 953–964 (2002)
- Sonnleitner E, Moll I, and Blasi U, Functional replacement of the *Escherichia coli* *hfq* gene by the homologue of *Pseudomonas aeruginosa*, *Microbiology* 148, 883–891 (2002)
- Soutourina O, Kolb A, Krin E, Laurent-Winter C, Rimsky S, Danchin A, and Bertin P, Multiple control of flagellum biosynthesis in *Escherichia coli*: role of H-NS protein and the cyclic AMP-catabolite activator protein complex in transcription of the *flhDC* master operon, *J Bacteriol* 181, 7500–7508 (1999)

- Soutourina OA, Semenova EA, Parfenova VV, Danchin A, and Bertin P, Control of bacterial motility by environmental factors in polarly flagellated and peritrichous bacteria isolated from Lake Baikal, *Appl Environ Microbiol* 67, 3852–3859 (2001)
- Srinivas PR, Srivastava S, Hanash S, and Wright GL Jr, Proteomics in early detection of cancer, *Clin Chem* 47, 1901–1911 (2001)
- Stanislaus R, Jiang LH, Swartz M, Arthur J, and Almeida JS, An XML standard for the dissemination of 2D gel electrophoresis data complemented with mass spectrometry results, *BMC Bioinformatics* 5, 9 (2004)
- Sturm R, Baumruker T, Franza BR Jr, and Herr W, A 100-kD HeLa cell octamer binding protein (OBP100) interacts differently with two separate octamer-related sequences within the SV40 enhancer, *Genes Dev* 1, 1147–1160 (1987)
- Sumpio BE, Banes AJ, Buckley M, and Johnson G Jr, Alterations in aortic endothelial cell morphology and cytoskeletal protein synthesis during cyclic tensile deformation, *J Vasc Surg* 7, 130–138 (1988)
- Sun J and Barbieri JT, *Pseudomonas aeruginosa* ExoT ADP-ribosylates CT10 regulator of kinase (Crk) proteins, *J Biol Chem* 278, 32794–32800 (2003)
- Sweatt JD and Kandel ER, Persistent and transcriptionally-dependent increase in protein phosphorylation in long-term facilitation of *Aplysia* sensory neurons, *Nature* 339, 51–54 (1989)
- Sweatt JD, Kennedy TE, Wager-Smith K, Gawinowicz MA, Barzilai A, Karl KA, and Kandel ER, Development of a database of amino acid sequences for proteins identified and isolated on two-dimensional polyacrylamide gels, *Electrophoresis* 10, 152–157 (1989)
- Takashima M, Kuramitsu Y, Yokoyama Y, Iizuka N, Toda T, Sakaida I, Okita K, Oka M, and Nakamura K, Proteomic profiling of heat shock protein 70 family members as biomarkers for hepatitis C virus-related hepatocellular carcinoma, *Proteomics* 3, 2487–2493 (2003)
- Taurin S, Seyrantepe V, Orlov SN, Tremblay TL, Thibault P, Bennett MR, Hamet P, and Pshzhetsky AV, Proteome analysis and functional expression identify mortalin as an antiapoptotic gene induced by elevation of $[Na^+]_i/[K^+]_i$ ratio in cultured vascular smooth muscle cells, *Circ Res* 91, 915–922 (2002)
- Taylor SW, Warnock DE, Glenn GM, Zhang B, Fahy E, Gaucher SP, Capaldi RA, Gibson BW, and Ghosh SS, An alternative strategy to determine the mitochondrial proteome using sucrose gradient fractionation and 1D PAGE on highly purified human heart mitochondria, *J Proteome Res* 1, 451–458 (2002)
- Terry DE and Desiderio DM, Between-gel reproducibility of the human cerebrospinal fluid proteome, *Proteomics* 3, 1962–1979 (2003)
- Toda T, Proteome and proteomics for the research on protein alterations in aging, *Ann NY Acad Sci* 928, 71–78 (2001)
- Toda T and Sugimoto M, Proteome analysis of Epstein-Barr virus-transformed B-lymphoblasts and the proteome database, *J Chromatogr B Analyt Technol Biomed Life Sci* 787, 197–206 (2003)
- Tryoen-Toth P, Richert S, Sohm B, Mine M, Marsac C, Van Dorsselaer A, Leize E, and Florentz C, Proteomic consequences of a human mitochondrial tRNA mutation beyond the frame of mitochondrial translation, *J Biol Chem* 278, 24314–24323 (2003)
- Turck N, Richert S, Gendry P, Stutzmann J, Kedinger M, Leize E, Simon-Assmann P, Van Dorsselaer A, and Launay JF, Proteomic analysis of nuclear proteins from proliferative and differentiated human colonic intestinal epithelial cells, *Proteomics* 4, 93–105 (2004)
- Tuttle JL, Hahn TL, Sanders BM, Witzmann FA, Miller SJ, Dalsing MC, and Unthank JL, Impaired collateral development in mature rats, *Am J Physiol Heart Circ Physiol* 283, H146–H155 (2002)
- ul Haq R and Chytil F, Early effects of retinol and retinoic acid on protein synthesis in retinol deficient rat testes, *Biochem Biophys Res Commun* 151, 53–60 (1988)
- Umar A, Ooms MP, Luider TM, Grootegoed JA, and Brinkmann AO, Proteomic profiling of epididymis and vas deferens: Identification of proteins regulated during rat genital tract development, *Endocrinology* 144, 4637–4647 (2003)
- Valadi H, Valadi A, Ansell R, Gustafsson L, Adler L, Norbeck J, and Blomberg A, NADH-reductive stress in *Saccharomyces cerevisiae* induces the expression of the minor isoform of glyceraldehyde-3-phosphate dehydrogenase (TDH1), *Curr Genet* 45, 90–95 (2004)
- van Anken E, Romijn EP, Maggioni C, Mezghrani A, Sitia R, Braakman I, and Heck AJ, Sequential waves of functionally related proteins are expressed when B cells prepare for antibody secretion, *Immunity* 18, 243–253 (2003)
- van der Vlies D, Pap EH, Post JA, Celis JE, and Wirtz KW, Endoplasmic reticulum resident proteins of normal human dermal fibroblasts are the major targets for oxidative stress induced by hydrogen peroxide, *Biochem J* 366, 825–830 (2002)
- Vanrobaeys F, Devreese B, Lecocq E, Rychlewski L, De Smet L, and Van Beeumen J, Proteomics of the dissimilatory iron-reducing bacterium *Shewanella oneidensis* MR-1, using a matrix-assisted laser desorption/ionization-tandem-time of flight mass spectrometer, *Proteomics* 3, 2249–2257 (2003)
- Vargas Mdel C, Encarnacion S, Davalos A, Reyes-Perez A, Mora Y, Garcia-de los Santos A, Brom S, and Mora J, Only one catalase, katG, is detectable in *Rhizobium etli*, and is encoded along with the regulator OxyR on a plasmid replicon, *Microbiology* 149, 1165–1176 (2003)
- Vecerek B, Moll I, Afonyushkin T, Kaberdin V, and Blasi U, Interaction of the RNA chaperone Hfq with mRNAs: direct and indirect roles of Hfq in iron metabolism of *Escherichia coli*, *Mol Microbiol* 50, 897–909 (2003)

- Vilain S, Cosette P, Zimmerlin I, Dupont JP, Junter GA, and Jouenne T, Biofilm proteome: homogeneity or versatility? *J Proteome Res* 3, 132–136 (2004)
- Vincek V, Nassiri M, Nadji M, and Morales AR, A tissue fixative that protects macromolecules (DNA, RNA, and protein) and histomorphology in clinical samples, *Lab Invest* 83, 1427–1435 (2003)
- Vivanco F, Lopez-Bescos L, Tunon J, and Egido J, Proteomics and cardiovascular disease, *Rev Esp Cardiol* 56, 289–302 (2003)
- Vohradsky J, Janda I, Grunenfelder B, Berndt P, Roder D, Langen H, Weiser J, and Jenal U, Proteome of *Caulobacter crescentus* cell cycle publicly accessible on SWICZ server, *Proteomics* 3, 1874–1882 (2003)
- Vohradsky J, Li XM, Dale G, Folcher M, Nguyen L, Viollier PH, and Thompson CJ, Developmental control of stress stimulons in *Streptomyces coelicolor* revealed by statistical analyses of global gene expression patterns, *J Bacteriol* 182, 4979–4986 (2000)
- Wagner WH, Henderson RM, Hicks HE, Banes AJ, and Johnson G Jr, Differences in morphology, growth rate, and protein synthesis between cultured arterial and venous endothelial cells, *J Vasc Surg* 8, 509–519 (1988)
- Wang W, Vignani R, Scali M, Sensi E, and Cresti M, Post-translational modifications of alpha-tubulin in *Zea mays* L. are highly tissue specific, *Planta* 218, 460–465 (2004)
- Watanabe H, Hamada H, Yamada N, Sohda S, Yamakawa-Kobayashi K, Yoshikawa H, and Arinami T, Proteome analysis reveals elevated serum levels of clusterin in patients with preeclampsia, *Proteomics* 4, 537–543 (2004)
- Wattiez R, Noel-Georis I, Cruyt C, Broeckeaert F, Bernard A, and Falmagne P, Susceptibility to oxidative stress: proteomic analysis of bronchoalveolar lavage from ozone-sensitive and ozone-resistant strains of mice, *Proteomics* 3, 658–665 (2003)
- Welch WJ, Garrels JI, Thomas GP, Lin JJ, and Feramisco JR, Biochemical characterization of the mammalian stress proteins and identification of two stress proteins as glucose- and Ca²⁺-ionophore-regulated proteins, *J Biol Chem* 258, 7102–7111 (1983)
- Wemekamp-Kamphuis HH, Karatzas AK, Wouters JA, and Abee T, Enhanced levels of cold shock proteins in *Listeria monocytogenes* LO28 upon exposure to low temperature and high hydrostatic pressure, *Appl Environ Microbiol* 68, 456–463 (2002)
- West KA, Yan L, Shadrach K, Sun J, Hasan A, Miyagi M, Crabb JS, Hollyfield JG, Marmorstein AD, and Crabb JW, Protein database, human retinal pigment epithelium, *Mol Cell Proteomics* 2, 37–49 (2003)
- Wheelock AM, Zhang L, Tran MU, Morin D, Penn S, Buckpitt AR, and Plopper CG, Isolation of rodent airway epithelial cell proteins facilitates in vivo proteomics studies of lung toxicity, *Am J Physiol Lung Cell Mol Physiol* 286, L399–L410 (2004)
- White MY, Cordwell SJ, McCarron HC, Tchen AS, Hambly BD, and Jeremy R, Modifications of myosin-regulatory light chain correlate with function of stunned myocardium, *J Mol Cell Cardiol* 35, 833–840 (2003)
- Wilke A, Ruckert C, Bartels D, Dondrup M, Goesmann A, Huser AT, Kespohl S, Linke B, Mahne M, McHardy A, Puhler A, and Meyer F, Bioinformatics support for high-throughput proteomics, *J Biotechnol* 106, 147–156 (2003)
- Wiseman RW, Lambert ME, Lamb PW, Garrels JI, and Barrett JC, Alterations in gene expression at various stages of neoplastic transformation of syrian hamster embryo (SHE) cells, in Moses HL, Lengyel P, Charles D and Stiles CD (eds) *Growth Inhibitory and Cytotoxic Polypeptides. Proceedings of a Genentech–Smith, Kline & French–Triton Biosciences–UCLA Symposium*, Keystone, Colorado, January 24–30, 1988, AR Liss, New York (1989)
- Witzmann FA and Li J, Cutting-edge technology. II. Proteomics: core technologies and applications in physiology, *Am J Physiol Gastrointest Liver Physiol* 282, G735–G741 (2002)
- Witzmann FA, Bobb A, Briggs GB, Coppage HN, Hess RA, Li J, Pedrick NM, Ritchie GD, Iii JR, and Still KR, Analysis of rat testicular protein expression following 91-day exposure to JP-8 jet fuel vapor, *Proteomics* 3, 1016–1027 (2003)
- Witzmann FA, Clack JW, Geiss K, Hussain S, Juhl MJ, Rice CM, and Wang C, Proteomic evaluation of cell preparation methods in primary hepatocyte cell culture, *Electrophoresis* 23, 2223–2232 (2002)
- Witzmann FA, Li J, Strother WN, McBride WJ, Hunter L, Crabb DW, Lumeng L, and Li TK, Innate differences in protein expression in the nucleus accumbens and hippocampus of inbred alcohol-preferring and nonpreferring rats, *Proteomics* 3, 1335–1344 (2003)
- Wu Y, Mou Z, Li J, Zhou W, Wei B, and Zou L, Identification of a S100 calcium-binding protein expressed in HL-60 cells treated with all-trans retinoic acid by two-dimensional electrophoresis and mass spectrometry, *Leuk Res* 28, 203–207 (2004)
- Xiao GG, Wang M, Li N, Loo JA, and Nel AE, Use of proteomics to demonstrate a hierarchical oxidative stress response to diesel exhaust particle chemicals in a macrophage cell line, *J Biol Chem* 278, 50781–50790 (2003)
- Xiong XD, Li EM, Xu LY, Chen HB, Chen L, Cai WJ, Han YL, Shen ZY, and Zeng Y, Separation and identification of differentially expressed nuclear matrix proteins between human esophageal immortalized and carcinomatous cell lines, *World J Gastroenterol* 9, 2143–2148 (2003)
- Xiong XD, Xu LY, Shen ZY, Cai WJ, Luo JM, Han YL, and Li EM, Identification of differentially expressed proteins between human esophageal immortalized and carcinomatous cell lines by two-dimensional electrophoresis and MALDI-TOF-mass spectrometry, *World J Gastroenterol* 8, 777–781 (2002)
- Yoon H, Lim S, Heu S, Choi S, and Ryu S, Proteome analysis of *Salmonella enterica* serovar Typhimurium *fts* mutant, *FEMS Microbiol Lett* 226, 391–396 (2003)

- Yu JH, Yun SY, Lim JW, Kim H, and Kim KH, Mass spectrometry and tandem mass spectrometry analysis of rat mitochondrial ATP synthase: up-regulation in pancreatic acinar cells treated with cerulein, *Proteomics* 3, 2437–2445 (2003)
- Yu JH, Yun SY, Lim JW, Kim H, and Kim KH, Proteome analysis of rat pancreatic acinar cells: implication for cerulein-induced acute pancreatitis, *Proteomics* 3, 2446–2453 (2003)
- Yu YL, Huang ZY, Yang PY, Rui YC, and Yang PY, Proteomic studies of macrophage-derived foam cell from human U937 cell line using two-dimensional gel electrophoresis and tandem mass spectrometry, *J Cardiovasc Pharmacol* 42, 782–789 (2003)
- Yu YL, Yang PY, Fan HZ, Huang ZY, Rui YC, and Yang PY, Protein expressions in macrophage-derived foam cells: comparative analysis by two-dimensional gel electrophoresis, *Acta Pharmacol Sin* 24, 873–877 (2003)
- Yuan X and Desiderio DM, Proteomics analysis of phosphotyrosyl-proteins in human lumbar cerebrospinal fluid, *J Proteome Res* 2, 476–487 (2003)
- Zhan X and Desiderio D, Heterogeneity analysis of the human pituitary proteome, *Clin Chem* 49, 1740–1751 (2003)
- Zhan X and Desiderio DM, Spot volume vs. amount of protein loaded onto a gel: A detailed, statistical comparison of two gel electrophoresis systems, *Electrophoresis* 24, 1818–1833 (2003)
- Zhan X and Desiderio D, Differences in the spatial and quantitative reproducibility between two second-dimensional gel electrophoresis systems, *Electrophoresis* 24, 1834–1846 (2003)
- Zhan X and Desiderio D, A reference map of a human pituitary adenoma proteome, *Proteomics* 3, 699–713 (2003)
- Zhan XQ, Guan YJ, Li C, Chen ZC, Xie JY, Chen P, and Liang SP, [Differential proteomic analysis of human lung adenocarcinoma cell line A-549 and of normal cell line HBE], *Sheng Wu Hua Xue Yu Sheng Wu Wu Li Xue Bao (Shanghai)* 34, 50–56 (2002)
- Zhang B, Su YP, Ai GP, Liu XH, Wang FC, and Cheng TM, Differentially expressed proteins of gamma-ray irradiated mouse intestinal epithelial cells by two-dimensional electrophoresis and MALDI-TOF mass spectrometry, *World J Gastroenterol* 9, 2726–2731 (2003)