

Odyssey® Publications

Selected from 251 publications in April 1, 2011 – June 30, 2011 on HighWire

Summer 2011

Volume 12

TABLE OF CONTENTS

	Page
Apoptosis	2
Autophagy	2
Angiogenesis	2
Cancer	3
Cardiovascular	4
Cell Adhesion, Migration and Invasion	5
Cell Cycle/DNA Replication	6
Cell Signaling	7
Diabetes/Metabolism/Obesity	8
Hypoxia	9
Immunology	9
Ion Channels	11
MicroRNA	11
Muscle	12
Neuroscience	12
Nephrology/Kidney	14
Phosphatases	14
Receptors	15
RNAi	16
Transcription Factors	17
Ubiquitination and Sumoylation	18
Virology/Microbiology	19
Applications	18-20
<i>In-Cell Western™ Assays</i>	
<i>IRDye® 800CW Toxicity Study</i>	
<i>Membrane Stripping</i>	
<i>Micro-Western Assay (Alternative to Reverse Phase Arrays)</i>	
<i>MousePOD® Accessories</i>	
<i>Reverse Phase Protein Arrays</i>	
<i>RNA/Protein Interactions</i>	
<i>siRNA Screening</i>	
<i>Western (Quantitative, Molecules/Cell)</i>	
PNAS and Science Publications This Quarter	23

LI-COR®

Experience Excellence

PUBLICATIONS REFERENCING ODYSSEY DATA

RESEARCH AREAS

APOPTOSIS

Hypersensitivity of mtDNA-depleted cells to staurosporine-induced apoptosis: roles of Bcl-2 downregulation and cathepsin B

Guillaume Rommelaere, Sébastien Michel, Ludovic Mercy, Antoine Fattaccioli, Catherine Demazy, Noelle Ninane, Andrée Houbion, Patricia Renard, and Thierry Arnould
Am J Physiol Cell Physiol, May 2011; 300: C1090 - C1106.

Stat3 signaling in acute myeloid leukemia: ligand-dependent and -independent activation and induction of apoptosis by a novel small-molecule Stat3 inhibitor

Michele S. Redell, Marcos J. Ruiz, Todd A. Alonzo, Robert B. Gerbing, and David J. Tweardy
Blood, May 2011; 117: 5701 - 5709.

Annexin A5-Conjugated Polymeric Micelles for Dual SPECT and Optical Detection of Apoptosis

Rui Zhang, Wei Lu, Xiaoxia Wen, Miao Huang, Min Zhou, Dong Liang, and Chun Li
J. Nucl. Med., Jun 2011; 52: 958 - 964.

Urocortin 1 modulates immunosignaling in a rat model of colitis via corticotropin-releasing factor receptor 2

Jen Chang, Melanie R. Adams, Matthew S. Clifton, Min Liao, Julia H. Brooks, Burcu Hasdemir, and Aditi Bhargava
Am J Physiol Gastrointest Liver Physiol, May 2011; 300: G884 - G894.

3'UTR-truncated *Hmga2* cDNA causes MPN-like hematopoiesis by conferring a clonal growth advantage at the level of HSC in mice

Kazuhiko Ikeda, Philip J. Mason, and Monica Bessler
Blood, Jun 2011; 117: 5860 - 5869.

BAX supports the mitochondrial network, promoting bioenergetics in nonapoptotic cells

Rebecca J. Boohaker, Ge Zhang, Adina Loosley Carlson, Kathleen N. Nemecek, and Annette R. Khaled
Am J Physiol Cell Physiol, Jun 2011; 300: C1466 - C1478.

AUTOPHAGY

Temporal Regulation of Intracellular Organelle Homeostasis in T Lymphocytes by Autophagy

Wei Jia and You-Wen He
J. Immunol., May 2011; 186: 5313 - 5322.

Rapamycin Reverses Cellular Phenotypes and Enhances Mutant Protein Clearance in Hutchinson-Gilford Progeria Syndrome Cells

Kan Cao, John J. Graziotto, Cecilia D. Blair, Joseph R. Mazzulli, Michael R. Erdos, Dimitri Krainc, and Francis S. Collins
Science Translational Medicine, Jun 2011; 3: 89ra58.

ANGIOGENESIS

AMP-activated protein kinase rescues the angiogenic functions of endothelial progenitor cells via manganese superoxide dismutase induction in type 1 diabetes

Xiao-Rong Wang, Ming-Wei Zhang, Dan-Dan Chen, Yun Zhang, and Alex F. Chen
Am J Physiol Endocrinol Metab, Jun 2011; 300: E1135 - E1145.

A Critical Role of Gβγ in Tumorigenesis and Metastasis of Breast Cancer

Xiaoyun Tang, Zhizeng Sun, Caitlin Runne, Joshua Madsen, Frederick Domann, Michael Henry, Fang Lin, and Songhai Chen
J. Biol. Chem., Apr 2011; 286: 13244 - 13254.

C/EBPβ Mediates Growth Hormone-Regulated Expression of Multiple Target Genes

Tracy X. Cui, Grace Lin, Christopher R. LaPensee, Anda-Alexandra Calinescu, Maanjot Rathore, Cale Streeter, Graciela Piwien-Pilipuk, Nathan Lanning, Hui Jin, Christin Carter-Su, Zhaohui S. Qin, and Jessica Schwartz
Mol. Endocrinol., Apr 2011; 25: 681 - 693.

Urocortin 1 modulates immunosignaling in a rat model of colitis via corticotropin-releasing factor receptor 2

Jen Chang, Melanie R. Adams, Matthew S. Clifton, Min Liao, Julia H. Brooks, Burcu Hasdemir, and Aditi Bhargava
Am J Physiol Gastrointest Liver Physiol, May 2011; 300: G884 - G894.

CANCER**Targeted Therapy of VEGFR2 and EGFR Significantly Inhibits Growth of Anaplastic Thyroid Cancer in an Orthotopic Murine Model**

Maria K. Gule, Yunyun Chen, Daisuke Sano, Mitchell J. Frederick, Ge Zhou, Mei Zhao, Zvonimir L. Milas, Chad E. Galer, Ying C. Henderson, Samar A. Jasser, David L. Schwartz, James A. Bankson, Jeffrey N. Myers, and Stephen Y. Lai
Clin. Cancer Res., Apr 2011; 17: 2281 - 2291.

Vandetanib Restores Head and Neck Squamous Cell Carcinoma Cells' Sensitivity to Cisplatin and Radiation *in vivo* and *in vitro*

Daisuke Sano, Fumihiko Matsumoto, David R. Valdecanas, Mei Zhao, David P. Molkenkine, Yoko Takahashi, Ehab Y. Hanna, Vali Papadimitrakopoulou, John Heymach, Luka Milas, and Jeffrey N. Myers
Clin. Cancer Res., Apr 2011; 17: 1815 – 1827.

High Plasma Levels and Effective Lymphatic Uptake of Docetaxel in an Orally Available Nanotransporter Formulation

Taher Nassar, Suha Attili-Qadri, Oshrat Harush-Frenkel, Shimon Farber, Shimon Lecht, Philip Lazarovici, and Simon Benita
Cancer Res., Apr 2011; 71: 3018 - 3028.

A Critical Role of G β in Tumorigenesis and Metastasis of Breast Cancer

Xiaoyun Tang, Zhizeng Sun, Caitlin Runne, Joshua Madsen, Frederick Domann, Michael Henry, Fang Lin, and Songhai Chen
J. Biol. Chem., Apr 2011; 286: 13244 - 13254.

A Novel HSP90 Inhibitor Delays Castrate-Resistant Prostate Cancer without Altering Serum PSA Levels and Inhibits Osteoclastogenesis

Francois Lamoureux, Christian Thomas, Min-Jean Yin, Hidetoshi Kuruma, Ladan Fazli, Martin E. Gleave, and Amina Zoubeidi
Clin. Cancer Res., Apr 2011; 17: 2301 - 2313.

Stromal LRP1 in Lung Adenocarcinoma Predicts Clinical Outcome

He Meng, Guoan Chen, Xiaojie Zhang, Zhuwen Wang, Dafydd G. Thomas, Thomas J. Giordano, David G. Beer, and Michael M. Wang
Clin. Cancer Res., Apr 2011; 17: 2426 - 2433.

High-Throughput Screen Identifies Novel Inhibitors of Cancer Biomarker α -Methylacyl Coenzyme A Racemase (AMACR/P504S)

Brice A.P. Wilson, Haofan Wang, Benjamin A. Nacev, Ronnie C. Mease, Jun O. Liu, Martin G. Pomper, and William B. Isaacs
Mol. Cancer Ther., May 2011; 10: 825 - 838.

Two-tiered Approach Identifies a Network of Cancer and Liver Disease-related Genes Regulated by miR-122

Daniel R. Boutz, Patrick J. Collins, Uthra Suresh, Mingzhu Lu, Cristina M. Ramírez, Carlos Fernández-Hernando, Yufei Huang, Raquel de Sousa Abreu, Shu-Yun Le, Bruce A. Shapiro, Angela M. Liu, John M. Luk, Shelley Force Aldred, Nathan D. Trinklein, Edward M. Marcotte, and Luiz O. F. Penalva
J. Biol. Chem., May 2011; 286: 18066 - 18078.

Phosphoglucose Isomerase/Autocrine Motility Factor Mediates Epithelial-Mesenchymal Transition Regulated by miR-200 in Breast Cancer Cells

Aamir Ahmad, Amro Aboukameel, Dejuan Kong, Zhiwei Wang, Seema Sethi, Wei Chen, Fazlul H. Sarkar, and Avraham Raz
Cancer Res., May 2011; 71: 3400 - 3409.

Pharmacological inhibition of sphingosine kinase isoforms alters estrogen receptor signaling in human breast cancer

James W Antoon, William D Meacham, Melyssa R Bratton, Evelyn M Slaughter, Lyndsay V Rhodes, Hasina B Ashe, Thomas E Wiese, Matthew E Burow, and Barbara S Beckman
J. Mol. Endocrinol., Jun 2011; 46: 205 - 216.

Mutationally Activated BRAF^{V600E} Elicits Papillary Thyroid Cancer in the Adult Mouse

Roch-Philippe Charles, Gioia Iezza, Elena Amendola, David Dankort, and Martin McMahon
Cancer Res., Jun 2011; 71: 3863 - 3871.

***In vivo* diagnosis of murine pancreatic intraepithelial neoplasia and early-stage pancreatic cancer by molecular imaging**

Stefan Eser, Marlena Messer, Philipp Eser, Alexander von Werder, Barbara Seidler, Monther Bajbouj, Roger Vogelmann, Alexander Meining, Johannes von Burstin, Hana Algül, Philipp Pagel, Angelika E. Schnieke, Irene Esposito, Roland M. Schmid, Günter Schneider, and Dieter Saur
PNAS, Jun 2011; 108: 9945 - 9950.

Epithelial Cell Organization Suppresses Myc Function by Attenuating Myc Expression

David R. Simpson, Min Yu, Siyuan Zheng, Zhongming Zhao, Senthil K. Muthuswamy, and William P. Tansey
Cancer Res., Jun 2011; 71: 3822 - 3830.

Antitumor and Antiangiogenic Activities of BMS-690514, an Inhibitor of Human EGF and VEGF Receptor Kinase Families

Tai W. Wong, Francis Y. Lee, Stuart Emanuel, Craig Fairchild, Joseph Fagnoli, Brian Fink, Ashvinikumar Gavai, Amy Hammell, Benjamin Henley, Christine Hilt, John T. Hunt, Bala Krishnan, Daniel Kukral, Anne Lewin, Harold Malone, Derek Norris, Simone Oppenheimer, Gregory Vite, and Chiang Yu
Clin. Cancer Res., Jun 2011; 17: 4031 - 4041.

HOXC9 Links Cell-Cycle Exit and Neuronal Differentiation and Is a Prognostic Marker in Neuroblastoma

Ling Mao, Jane Ding, Yunhong Zha, Liqun Yang, Brian A. McCarthy, William King, Hongjuan Cui, and Han-Fei Ding
Cancer Res., Jun 2011; 71: 4314 - 4324.

Novel Changes in Glycosylation of Serum Apo-J in Patients with Hepatocellular Carcinoma

Mary Ann Comunale, Mengjun Wang, Lucy Rodemich-Betesh, Julie Hafner, Anne Lamontagne, Andrew Klein, Jorge Marrero, Adrian M. Di Bisceglie, Robert Gish, Timothy Block, and Anand Mehta
Cancer Epidemiol. Biomarkers Prev., Jun 2011; 20: 1222 - 1229.

Mutationally Activated BRAF^{V600E} Elicits Papillary Thyroid Cancer in the Adult Mouse

Roch-Philippe Charles, Gioia Iezza, Elena Amendola, David Dankort, and Martin McMahon
Cancer Res., Jun 2011; 71: 3863 - 3871.

Cardiovascular

Tumor Necrosis Factor- α Accelerates the Calcification of Human Aortic Valve Interstitial Cells Obtained from Patients with Calcific Aortic Valve Stenosis via the BMP2-Dlx5 Pathway

Zaiqiang Yu, Kazuhiko Seya, Kazuyuki Daitoku, Shigeru Motomura, Ikuo Fukuda, and Ken-Ichi Furukawa
J. Pharmacol. Exp. Ther., Apr 2011; 337: 16 - 23.

Soluble Form of the (Pro)Renin Receptor Is Augmented in the Collecting Duct and Urine of Chronic Angiotensin II-Dependent Hypertensive Rats

Alexis A. Gonzalez, Lucienne S. Lara, Christina Luffman, Dale M. Seth, and Minolfa C. Prieto
Hypertension, Apr 2011; 57: 859 - 864.

Inhibition of renal caveolin-1 reduces natriuresis and produces hypertension in sodium-loaded rats

John J. Gildea, Brandon A. Kemp, Nancy L. Howell, Robert E. Van Sciver, Robert M. Carey, and Robin A. Felder
Am J Physiol Renal Physiol, Apr 2011; 300: F914 - F920.

Antisense oligonucleotide reduction of apoB-ameliorated atherosclerosis in LDL receptor-deficient mice

Adam E. Mullick, Wuxia Fu, Mark J. Graham, Richard G. Lee, Donna Witchell, Thomas A. Bell, Charles P. Whipple, and Rosanne M. Crooke
J. Lipid Res., May 2011; 52: 885 - 896.

Induction of Heme Oxygenase 1 Attenuates Placental Ischemia-Induced Hypertension

Eric M. George, Kathy Cockrell, Marietta Aranay, Eva Csongradi, David E. Stec, and Joey P. Granger
Hypertension, May 2011; 57: 941 - 948.

Changes in cross-bridge cycling underlie muscle weakness in patients with tropomyosin 3-based myopathy

Coen A.C. Ottenheijm, Michael W. Lawlor, Ger J.M. Stienen, Henk Granzier, and Alan H. Beggs

Hum. Mol. Genet., May 2011; 20: 2015 - 2025.**Myocardin is differentially required for the development of smooth muscle cells and cardiomyocytes**

Mark H. Hoofnagle, Ronald L. Nepl, Erica L. Berzin, G. C. Teg Pipes, Eric N. Olson, Brian W. Wamhoff, Avril V. Somlyo, and Gary K. Owens

Am J Physiol Heart Circ Physiol, May 2011; 300: H1707 - H1721.**Protein Kinase C β_{II} -Mediated Phosphorylation of Endothelial Nitric Oxide Synthase Threonine 495 Mediates the Endothelial Dysfunction Induced by FK506 (Tacrolimus)**

Valorie L. Chiasson, Matthew A. Quinn, Kristina J. Young, and Brett M. Mitchell

J. Pharmacol. Exp. Ther., Jun 2011; 337: 718 - 723.**Cardiac-specific mindin overexpression attenuates cardiac hypertrophy via blocking AKT/GSK3 β and TGF- β 1-Smad signalling**

Ling Yan, Xiang Wei, Qi-Zhu Tang, Jinghua Feng, Yan Zhang, Chen Liu, Zhou-Yan Bian, Lian-Feng Zhang, Manyin Chen, Xue Bai, Ai-Bing Wang, John Fassett, Yingjie Chen, You-Wen He, Qinglin Yang, Peter P. Liu, and Hongliang Li

Cardiovasc Res, Jun 2011; 10.1093/cvr/cvr159.**Phosphatase-Resistant Gap Junctions Inhibit Pathological Remodeling and Prevent Arrhythmias**

Benjamin F. Remo, Jiayang Qu, Frank M. Volpicelli, Steven Giovannone, Daniel Shin, Joshua Lader,

Fang-yu Liu, Jie Zhang, Danielle S. Lent, Gregory E. Morley, and Glenn I. Fishman

Circ. Res., Jun 2011; 108: 1459 - 1466.

Cell Adhesion, Migration and Invasion

Macrophage Migration Inhibitory Factor and CD74 Regulate Macrophage Chemotactic Responses via MAPK and Rho GTPase

Huapeng Fan, Pam Hall, Leilani L. Santos, Julia L. Gregory, Gunter Fingerle-Rowson, Richard Bucala, Eric F. Morand, and Michael J. Hickey

J. Immunol., Apr 2011; 186: 4915 - 4924.**T-bet is induced by interferon- γ to mediate chemokine secretion and migration in human airway smooth muscle cells**

Cherie A. Singer

Am J Physiol Lung Cell Mol Physiol, Apr 2011; 300: L633 - L641.**CD43 interaction with ezrin-radixin-moesin (ERM) proteins regulates T-cell trafficking and CD43 phosphorylation**

J. L. Cannon, P. D. Mody, K. M. Blaine, E. J. Chen, A. D. Nelson, L. J. Sayles, T. V. Moore, B. S. Clay, N. O. Dulin, R. A. Shilling, J. K. Burkhardt, and A. I. Sperling

Mol. Biol. Cell, Apr 2011; 22: 954 - 963.**Hematopoietic Lineage Cell-Specific Protein 1 Functions in Concert with the Wiskott-Aldrich Syndrome Protein To Promote Podosome Array Organization and Chemotaxis in Dendritic Cells**

Deborah A. Klos Dehring, Fiona Clarke, Brendon G. Ricart, Yanping Huang, Timothy S. Gomez, Edward K. Williamson, Daniel A. Hammer, Daniel D. Billadeau, Yair Argon, and Janis K. Burkhardt

J. Immunol., Apr 2011; 186: 4805 - 4818.**Mechanistic Control of Carcinoembryonic Antigen-related Cell Adhesion Molecule-1 (CEACAM1) Splice Isoforms by the Heterogeneous Nuclear Ribonuclear Proteins hnRNP L, hnRNP A1, and hnRNP M**

Kenneth J. Dery, Shikha Gaur, Marieta Gencheva, YunYen, John E. Shively, and Rajesh K. Gaur

J. Biol. Chem., May 2011; 286: 16039 - 16051.**Hax1 regulates neutrophil adhesion and motility through RhoA**

Peter J. Cavnar, Erwin Berthier, David J. Beebe, and Anna Huttenlocher

J. Cell Biol., May 2011; 193: 465 - 473.

Phosphoglucose Isomerase/Autocrine Motility Factor Mediates Epithelial-Mesenchymal Transition Regulated by miR-200 in Breast Cancer Cells

Aamir Ahmad, Amro Aboukameel, Dejuan Kong, Zhiwei Wang, Seema Sethi, Wei Chen, Fazlul H. Sarkar, and Avraham Raz
Cancer Res., May 2011; 71: 3400 - 3409.

Tropomyosin variants describe distinct functional subcellular domains in differentiated vascular smooth muscle cells

Cynthia Gallant, Sarah Appel, Philip Graceffa, Paul Leavis, Jim Jung-Ching Lin, Peter W. Gunning, Galina Schevzov, Christine Chaponnier, Jon DeGnore, William Lehman, and Kathleen G. Morgan
Am J Physiol Cell Physiol, Jun 2011; 300: C1356 - C1365.

Targeted Identification of Metastasis-associated Cell-surface Sialoglycoproteins in Prostate Cancer

Lifang Yang, Julius O. Nyalwidhe, Siqi Guo, Richard R. Drake, and O. John Semmes
Mol. Cell. Proteomics, Jun 2011; 10: M110.007294.

Regulation of membrane-cytoskeletal interactions by tyrosine phosphorylation of erythrocyte band 3

Emanuela Ferru, Katie Giger, Antonella Pantaleo, Estela Campanella, Jesse Grey, Ken Ritchie, Rosa Vono, Francesco Turrini, and Philip S. Low
Blood, Jun 2011; 117: 5998 - 6006.

CYLD Inhibits Tumorigenesis and Metastasis by Blocking JNK/AP1 Signaling at Multiple Levels

Paula Miliiani de Marval, Shazia Lutfeali, Jane Y. Jin, Benjamin Leshin, M. Angelica Selim, and Jennifer Y. Zhang
Cancer Prevention Research, Jun 2011; 4: 851 - 859.

***In vivo* functional analysis reveals specific roles for the integrin-binding sites of talin**

Stephanie J. Ellis, Mary Pines, Michael J. Fairchild, and Guy Tanentzapf
J. Cell Sci., Jun 2011; 124: 1844 - 1856.

Cell Cycle/DNA Replication

Human NDR Kinases Control G₁/S Cell Cycle Transition by Directly Regulating p21 Stability

Hauke Cornils, Reto S. Kohler, Alexander Hergovich, and Brian A. Hemmings
Mol. Cell. Biol., Apr 2011; 31: 1382 - 1395.

USP7/HAUSP stimulates repair of oxidative DNA lesions

Svetlana V. Khoronenkova, Irina I. Dianova, Jason L. Parsons, and Grigory L. Dianov
Nucleic Acids Res., Apr 2011; 39: 2604 - 2609.

Tumour suppressor ING1b maintains genomic stability upon replication stress

Ronald P. C. Wong, Hanyang Lin, Shahram Khosravi, Brad Piche, Seyed Mehdi Jafarnejad, David W. C. Chen, and Gang Li
Nucleic Acids Res., May 2011; 39: 3632 - 3642.

Activities of ICP0 Involved in the Reversal of Silencing of Quiescent Herpes Simplex Virus 1

Michael W. Ferenczy, Daniel J. Ranayhossaini, and Neal A. DeLuca
J. Virol., May 2011; 85: 4993 - 5002.

HOXC9 Links Cell-Cycle Exit and Neuronal Differentiation and Is a Prognostic Marker in Neuroblastoma

Ling Mao, Jane Ding, Yunhong Zha, Liqun Yang, Brian A. McCarthy, William King, Hongjuan Cui, and Han-Fei Ding
Cancer Res., Jun 2011; 71: 4314 - 4324.

Formation of a Stable RuvA Protein Double Tetramer Is Required for Efficient Branch Migration *in vitro* and for Replication Fork Reversal *in vivo*

Alison S. Bradley, Zeynep Baharoglu, Andrew Niewiarowski, Bénédicte Michel, and Irina R. Tsaneva
J. Biol. Chem., Jun 2011; 286: 22372 - 22383.

Direct regulation of Treslin by cyclin-dependent kinase is essential for the onset of DNA replication

Akiko Kumagai, Anna Shevchenko, Andrej Shevchenko, and William G. Dunphy
J. Cell Biol., Jun 2011; 193: 995 - 1007.

Impairment of Human Immunodeficiency Virus Type-1 Integrase SUMOylation Correlates with an Early Replication Defect

Alessia Zamborlini, Audrey Coiffic, Guillaume Beauclair, Olivier Delelis, Joris Paris, Yashuiro Koh, Fabian Magne, Marie-Lou Giron, Joelle Tobaly-Tapiero, Eric Deprez, Stephane Emiliani, Alan Engelman, Hugues deThé, and Ali Saïb
J. Biol. Chem., Jun 2011; 286: 21013 - 21022.

Interleukin-13 inhibits proliferation and enhances contractility of human airway smooth muscle cells without change in contractile phenotype

Paul-André Risse, Taisuke Jo, Fernando Suarez, Nobuaki Hirota, Barbara Tolloczko, Pasquale Ferraro, Peter Grutter, and James G. Martin
Am J Physiol Lung Cell Mol Physiol, Jun 2011; 300: L958 - L966.

Cell Signaling

A Critical Role of Gβγ in Tumorigenesis and Metastasis of Breast Cancer

Xiaoyun Tang, Zhizeng Sun, Caitlin Runne, Joshua Madsen, Frederick Domann, Michael Henry, Fang Lin, and Songhai Chen
J. Biol. Chem., Apr 2011; 286: 13244 - 13254.

Upregulation of ABCG2 by Romidepsin via the Aryl Hydrocarbon Receptor Pathway

Kenneth K.W. To, Robert Robey, Zhirong Zhan, Lois Bangiolo, and Susan E. Bates
Mol. Cancer Res., Apr 2011; 9: 516 - 527.

Notch1-mediated Signaling Induces MHC Class II Expression through Activation of Class II Transactivator Promoter III in Mast Cells

Nobuhiro Nakano, Chiharu Nishiyama, Hideo Yagita, Akemi Koyanagi, Hideoki Ogawa, and Ko Okumura
J. Biol. Chem., Apr 2011; 286: 12042 - 12048.

Lysate Microarrays Enable High-throughput, Quantitative Investigations of Cellular Signaling

Mark Sevecka, Alejandro Wolf-Yadlin, and Gavin MacBeath
Mol. Cell. Proteomics, Apr 2011; 10: M110.005363.

Upregulation of ABCG2 by Romidepsin via the Aryl Hydrocarbon Receptor Pathway

Kenneth K.W. To, Robert Robey, Zhirong Zhan, Lois Bangiolo, and Susan E. Bates
Mol. Cancer Res., Apr 2011; 9: 516 - 527.

STAT3 Mediates Resistance to MEK Inhibitor through MicroRNA miR-17

Bingbing Dai, Jieru Meng, Michael Peyton, Luc Girard, William G. Bornmann, Lin Ji, John D. Minna, Bingliang Fang, and Jack A. Roth
Cancer Res., May 2011; 71: 3658 - 3668.

Stat3 signaling in acute myeloid leukemia: ligand-dependent and -independent activation and induction of apoptosis by a novel small-molecule Stat3 inhibitor

Michele S. Redell, Marcos J. Ruiz, Todd A. Alonzo, Robert B. Gerbing, and David J. Tweardy
Blood, May 2011; 117: 5701 - 5709.

A Peptide-Based Target Screen Implicates the Protein Kinase CK2 in the Global Regulation of Caspase Signaling

James S. Duncan, Jacob P. Turowec, Kelly E. Duncan, Greg Vilks, Chenggang Wu, Bernhard Lüscher, Shawn S.-C. Li, Greg B. Gloor, and David W. Litchfield
Sci. Signal., May 2011; 4: ra30.

Acidic nuclear phosphoprotein 32kDa (ANP32)B-deficient mouse reveals a hierarchy of ANP32 importance in mammalian development

Patrick T. Reilly, Samia Afzal, Chiara Gorrini, Koren Lui, Yury V. Bukhman, Andrew Wakeham, Jillian Haight, Teo Wei Ling, Carol C. Cheung, Andrew J. Elia, Patricia V. Turner, and Tak Wah Mak
PNAS, Jun 2011; 108: 10243 - 10248.

Pharmacological inhibition of sphingosine kinase isoforms alters estrogen receptor signaling in human breast cancer

James W Antoon, William D Meacham, Melyssa R Bratton, Evelyn M Slaughter, Lyndsay V Rhodes, Hasina B Ashe, Thomas E Wiese, Matthew E Burow, and Barbara S Beckman
J. Mol. Endocrinol., Jun 2011; 46: 205 - 216.

Regulation of mammalian Gli proteins by Costal 2 and PKA in Drosophila reveals Hedgehog pathway conservation

Steven A. Marks and Daniel Kalderon
Development, Jun 2011; 138: 2533 - 2542.

CYLD Inhibits Tumorigenesis and Metastasis by Blocking JNK/AP1 Signaling at Multiple Levels

Paula Miliiani de Marval, Shazia Lutfeali, Jane Y. Jin, Benjamin Leshin, M. Angelica Selim, and Jennifer Y. Zhang
Cancer Prevention Research, Jun 2011; 4: 851 - 859.

Pharmacological inhibition of sphingosine kinase isoforms alters estrogen receptor signaling in human breast cancer

James W Antoon, William D Meacham, Melyssa R Bratton, Evelyn M Slaughter, Lyndsay V Rhodes, Hasina B Ashe, Thomas E Wiese, Matthew E Burow, and Barbara S Beckman
J. Mol. Endocrinol., Jun 2011; 46: 205 - 216.

Regulation of Src Family Kinases Involved in T Cell Receptor Signaling by Protein-tyrosine Phosphatase CD148

Ondrej Stepanek, Tomas Kalina, Peter Draber, Tereza Skopcova, Karel Svojgr, Pavla Angelisova, Vaclav Horejsi, Arthur Weiss, and Tomas Brdicka
J. Biol. Chem., Jun 2011; 286: 22101 - 22112.

Diabetes/ Metabolism/Obesity

Central Endocannabinoid Signaling Regulates Hepatic Glucose Production and Systemic Lipolysis

James D. O'Hare, Elizabeth Zielinski, Bob Cheng, Thomas Scherer, and Christoph Buettner
Diabetes, Apr 2011; 60: 1055 - 1062.

MicroRNA-29c Is a Signature MicroRNA under High Glucose Conditions That Targets Sprouty Homolog 1, and Its *in Vivo* Knockdown Prevents Progression of Diabetic Nephropathy

Jianyin Long, Yin Wang, Wenjian Wang, Benny H. J. Chang, and Farhad R. Danesh
J. Biol. Chem., Apr 2011; 286: 11837 - 11848.

Depletion of PINK1 affects mitochondrial metabolism, calcium homeostasis and energy maintenance

Bavo Heeman, Chris Van den Haute, Sarah-Ann Aelvoet, Federica Valsecchi, Richard J. Rodenburg, Veerle Reumers, Zeger Debyser, Geert Callewaert, Werner J. H. Koopman, Peter H. G. M. Willems, and Veerle Baekelandt
J. Cell Sci., Apr 2011; 124: 1115 - 1125.

Nutrient-Sensing Hypothalamic TXNIP Links Nutrient Excess to Energy Imbalance in Mice

Clémence Blouet and Gary J. Schwartz
J. Neurosci., Apr 2011; 31: 6019 - 6027.

Novel Fat Depot-Specific Mechanisms Underlie Resistance to Visceral Obesity and Inflammation in 11 β -Hydroxysteroid Dehydrogenase Type 1-Deficient Mice

Malgorzata Wamil, Jenny H. Battle, Sophie Turban, Tiina Kipari, David Seguret, Ricardo de Sousa Peixoto, Yvonne B. Nelson, Dominika Nowakowska, David Ferenbach, Lynne Ramage, Karen E. Chapman, Jeremy Hughes, Donald R. Dunbar, Jonathan R. Seckl, and Nicholas M. Morton
Diabetes, Apr 2011; 60: 1158 - 1167.

Role of Inflammation and Insulin Resistance in Endothelial Progenitor Cell Dysfunction

Cyrus V. Desouza, Frederick G. Hamel, Keshore Bidasee, and Kelly O'Connell
Diabetes, Apr 2011; 60: 1286 - 1294.

CD36 Protein Is Involved in Store-operated Calcium Flux, Phospholipase A2 Activation, and Production of Prostaglandin E2

Ondrej Kuda, Christopher M. Jenkins, James R. Skinner, Sung Ho Moon, Xiong Su, Richard W. Gross, and Nada A. Abumrad
J. Biol. Chem., May 2011; 286: 17785 - 17795.

Improved Insulin Sensitivity by GLUT12 Overexpression in Mice

Scott H. Purcell, Lauren B. Aerni-Flessner, Alexandra R. Willcockson, Kelly A. Diggs-Andrews, Simon J. Fisher, and Kelle H. Moley
Diabetes, May 2011; 60: 1478 - 1482.

Inhibition of Human Cytochrome P450 3A4 by Cholesterol

Raku Shinkyo and F. Peter Guengerich
J. Biol. Chem., May 2011; 286: 18426 - 18433.

CYP2C9-Mediated Metabolic Activation of Losartan Detected by a Highly Sensitive Cell-Based Screening Assay

Atsushi Iwamura, Tatsuki Fukami, Hiroko Hosomi, Miki Nakajima, and Tsuyoshi Yokoi
Drug Metab. Dispos., May 2011; 39: 838 - 846.

AMP-activated protein kinase rescues the angiogenic functions of endothelial progenitor cells via manganese superoxide dismutase induction in type 1 diabetes

Xiao-Rong Wang, Ming-Wei Zhang, Dan-Dan Chen, Yun Zhang, and Alex F. Chen
Am J Physiol Endocrinol Metab, Jun 2011; 300: E1135 - E1145.

Continuous 24-h nicotinic acid infusion in rats causes FFA rebound and insulin resistance by altering gene expression and basal lipolysis in adipose tissue

YoungTaek Oh, Ki-Sook Oh, Yong Min Choi, Anne Jokiah, Casey Donovan, Sangdun Choi, Insug Kang, and Jang H. Youn
Am J Physiol Endocrinol Metab, Jun 2011; 300: E1012 - E1021.

Separation of Fast from Slow Anabolism by Site-specific PEGylation of Insulin-like Growth Factor I (IGF-I)

Friedrich Metzger, Waseem Sajid, Stefanie Saenger, Christian Staudenmaier, Chris van der Poel, Bettina Sobottka, Angelika Schuler, Mandy Sawitzky, Raphael Poirier, Dietrich Tuerck, Eginhard Schick, Andreas Schaubmar, Friederike Hesse, Kurt Amrein, Hansruedi Loetscher, Gordon S. Lynch, Andreas Hoeflich, Pierre De Meyts, and Hans-Joachim Schoenfeld
J. Biol. Chem., Jun 2011; 286: 19501 - 19510.

A Potential Role for Angiotensin 2 in the Regulation of the Blood–Retinal Barrier in Diabetic Retinopathy

Sampathkumar Rangasamy, Ramprasad Srinivasan, Joann Maestas, Paul G. McGuire, and Arup Das
Invest. Ophthalmol. Vis. Sci., Jun 2011; 52: 3784 - 3791.

BAX supports the mitochondrial network, promoting bioenergetics in nonapoptotic cells

Rebecca J. Boohaker, Ge Zhang, Adina Loosley Carlson, Kathleen N. Nemecek, and Annette R. Khaled
Am J Physiol Cell Physiol, Jun 2011; 300: C1466 - C1478.

Hypoxia

Nicotinamide Mononucleotide Adenylyltransferase Is a Stress Response Protein Regulated by the Heat Shock Factor/Hypoxia-inducible Factor 1 α Pathway

Yousuf O. Ali, Ryan McCormack, Andrew Darr, and R. Grace Zhai
J. Biol. Chem., May 2011; 286: 19089 - 19099.

Mitochondrial c-Jun N-terminal Kinase (JNK) Signaling Initiates Physiological Changes Resulting in Amplification of Reactive Oxygen Species Generation

Jeremy W. Chambers and Philip V. LoGrasso
J. Biol. Chem., May 2011; 286: 16052 - 16062.

Immunology

Role of Inflammation and Insulin Resistance in Endothelial Progenitor Cell Dysfunction

Cyrus V. Desouza, Frederick G. Hamel, Keshore Bidasee, and Kelly O'Connell
Diabetes, Apr 2011; 60: 1286 - 1294.

Mal Mediates TLR-Induced Activation of CREB and Expression of IL-10

Mark Mellett, Paola Atzei, Ruaidhri Jackson, Luke A. O'Neill, and Paul N. Moynagh
J. Immunol., Apr 2011; 186: 4925 - 4935.

CD43 interaction with ezrin-radixin-moesin (ERM) proteins regulates T-cell trafficking and CD43 phosphorylation

J. L. Cannon, P. D. Mody, K. M. Blaine, E. J. Chen, A. D. Nelson, L. J. Sayles, T. V. Moore, B. S. Clay, N. O. Dulin, R. A. Shilling, J. K. Burkhardt, and A. I. Sperling
Mol. Biol. Cell, Apr 2011; 22: 954 - 963.

Modulation of Innate and Adaptive Immune Responses by Tofacitinib (CP-690,550)

Kamran Ghoreschi, Michael I. Jesson, Xiong Li, Jamie L. Lee, Sarbani Ghosh, Jason W. Alsup, James D. Warner, Masao Tanaka, Scott M. Steward-Tharp, Massimo Gadina, Craig J. Thomas, John C. Minnerly, Chad E. Storer, Timothy P. LaBranche, Zaher A. Radi, Martin E. Dowty, Richard D. Head, Debra M. Meyer, Nandini Kishore, and John J. O'Shea
J. Immunol., Apr 2011; 186: 4234 - 4243.

Hematopoietic Lineage Cell-Specific Protein 1 Functions in Concert with the Wiskott–Aldrich Syndrome Protein To Promote Podosome Array Organization and Chemotaxis in Dendritic Cells

Deborah A. Klos Dehring, Fiona Clarke, Brendon G. Ricart, Yanping Huang, Timothy S. Gomez, Edward K. Williamson, Daniel A. Hammer, Daniel D. Billadeau, Yair Argon, and Janis K. Burkhardt
J. Immunol., Apr 2011; 186: 4805 - 4818.

Macrophage Migration Inhibitory Factor and CD74 Regulate Macrophage Chemotactic Responses via MAPK and Rho GTPase

Huapeng Fan, Pam Hall, Leilani L. Santos, Julia L. Gregory, Gunter Fingerle-Rowson, Richard Bucala, Eric F. Morand, and Michael J. Hickey
J. Immunol., Apr 2011; 186: 4915 - 4924.

Complement System Dysregulation and Inflammation in the Retinal Pigment Epithelium of a Mouse Model for Stargardt Macular Degeneration

Roxana A. Radu, Jane Hu, Quan Yuan, Darcy L. Welch, Jacob Makshanoff, Marcia Lloyd, Stephen McMullen, Gabriel H. Travis, and Dean Bok
J. Biol. Chem., May 2011; 286: 18593 - 18601.

Urocortin 1 modulates immunosignaling in a rat model of colitis via corticotropin-releasing factor receptor 2

Jen Chang, Melanie R. Adams, Matthew S. Clifton, Min Liao, Julia H. Brooks, Burcu Hasdemir, and Aditi Bhargava
Am J Physiol Gastrointest Liver Physiol, May 2011; 300: G884 - G894.

The Pleckstrin Homology Domain in the SKAP55 Adapter Protein Defines the Ability of the Adapter Protein ADAP To Regulate Integrin Function and NF- κ B Activation

Brandon J. Burbach, Rupa Srivastava, Melissa A. Ingram, Jason S. Mitchell, and Yoji Shimizu
J. Immunol., Jun 2011; 186: 6227 - 6237.

p53 Negatively Regulates RGS13 Protein Expression in Immune Cells

Shoko Iwaki, Yunbiao Lu, Zhihui Xie, and Kirk M. Druey
J. Biol. Chem., Jun 2011; 286: 22219 - 22226.

Overexpression of the Dominant-Negative Form of Interferon Regulatory Factor 1 in Oligodendrocytes Protects against Experimental Autoimmune Encephalomyelitis

Zihua Ren, Yan Wang, Duan Tao, David Liebenson, Thomas Liggett, Rajendra Goswami, Robert Clarke, Dusan Stefoski, and Roumen Balabanov
J. Neurosci., Jun 2011; 31: 8329 - 8341.

Regulation of Src Family Kinases Involved in T Cell Receptor Signaling by Protein-tyrosine Phosphatase CD148

Ondrej Stepanek, Tomas Kalina, Peter Draber, Tereza Skopcova, Karel Svojgr, Pavla Angelisova, Vaclav Horejsi, Arthur Weiss, and Tomas Brdicka
J. Biol. Chem., Jun 2011; 286: 22101 - 22112.

Tight Control of STAT5 Activity Determines Human CD34-Derived Interstitial Dendritic Cell and Langerhans Cell Development

Lianne van de Laar, Aniek van den Bosch, Albertus T. J. Wierenga, Harry L. A. Janssen, Paul J. Coffey, and Andrea M. Woltman
J. Immunol., Jun 2011; 186: 7016 - 7024.

Ion Channels

Large-scale phosphoproteomic analysis of membrane proteins in renal proximal and distal tubule

Marina Feric, Boyang Zhao, Jason D. Hoffert, Trairak Pisitkun, and Mark A. Knepper
Am J Physiol Cell Physiol, Apr 2011; 300: C755 - C770.

Small GTPase Rab11b regulates degradation of surface membrane L-type Ca_v1.2 channels

Jabe M. Best, Jason D. Foell, Courtney R. Buss, Brian P. Delisle, Ravi C. Balijepalli, Craig T. January, and Timothy J. Kamp
Am J Physiol Cell Physiol, May 2011; 300: C1023 - C1033.

Rat diaphragm mitochondria have lower intrinsic respiratory rates than mitochondria in limb muscles

Mary L. Garcia-Cazarin, Jorge L. Gamboa, and Francisco H. Andrade
Am J Physiol Regulatory Integrative Comp Physiol, Jun 2011; 300: R1311 - R1315.

Cortactin Is Required for N-cadherin Regulation of Kv1.5 Channel Function

Lan Cheng, Aaron Yung, Manuel Covarrubias, and Glenn L. Radice
J. Biol. Chem., Jun 2011; 286: 20478 - 20489.

Phosphatase-Resistant Gap Junctions Inhibit Pathological Remodeling and Prevent Arrhythmias

Benjamin F. Remo, Jiaxiang Qu, Frank M. Volpicelli, Steven Giovannone, Daniel Shin, Joshua Lader, Fang-yu Liu, Jie Zhang, Danielle S. Lent, Gregory E. Morley, and Glenn I. Fishman
Circ. Res., Jun 2011; 108: 1459 - 1466.

MicroRNA

MicroRNA-29c Is a Signature MicroRNA under High Glucose Conditions That Targets Sprouty Homolog 1, and Its *in vivo* Knockdown Prevents Progression of Diabetic Nephropathy

Jianyin Long, Yin Wang, Wenjian Wang, Benny H. J. Chang, and Farhad R. Danesh
J. Biol. Chem., Apr 2011; 286: 11837 - 11848.

Hormonal Regulation of Sertoli Cell Micro-RNAs at Spermiation

Peter K. Nicholls, Craig A. Harrison, Kelly L. Walton, Robert I. McLachlan, Liza O'Donnell, and Peter G. Stanton
Endocrinology, Apr 2011; 152: 1670 - 1683.

Two-tiered Approach Identifies a Network of Cancer and Liver Disease-related Genes Regulated by miR-122

Daniel R. Boutz, Patrick J. Collins, Uthra Suresh, Mingzhu Lu, Cristina M. Ramirez, Carlos Fernández-Hernando, Yufei Huang, Raquel de Sousa Abreu, Shu-Yun Le, Bruce A. Shapiro, Angela M. Liu, John M. Luk, Shelley Force Aldred, Nathan D. Trinklein, Edward M. Marcotte, and Luiz O. F. Penalva
J. Biol. Chem., May 2011; 286: 18066 - 18078.

Coordinated Regulation of Extracellular Matrix Synthesis by the MicroRNA-29 Family in the Trabecular Meshwork

Guadalupe Villarreal, Jr, Dong-Jin Oh, Min Hyung Kang, and Douglas J. Rhee
Invest. Ophthalmol. Vis. Sci., May 2011; 52: 3391 - 3397.

STAT3 Mediates Resistance to MEK Inhibitor through MicroRNA miR-17

Bingbing Dai, Jieru Meng, Michael Peyton, Luc Girard, William G. Bornmann, Lin Ji, John D. Minna, Bingliang Fang, and Jack A. Roth
Cancer Res., May 2011; 71: 3658 - 3668.

Phosphoglucose Isomerase/Autocrine Motility Factor Mediates Epithelial-Mesenchymal Transition Regulated by miR-200 in Breast Cancer Cells

Aamir Ahmad, Amro Aboukameel, Dejuan Kong, Zhiwei Wang, Seema Sethi, Wei Chen, Fazlul H. Sarkar, and Avraham Raz
Cancer Res., May 2011; 71: 3400 - 3409.

Platelet microRNA-mRNA coexpression profiles correlate with platelet reactivity

Srikanth Nagalla, Chad Shaw, Xianguo Kong, Altaf A. Kondkar, Leonard C. Edelstein, Lin Ma, Junmei Chen, G. Stanley McKnight, José A. López, Linghai Yang, Ying Jin, Molly S. Bray, Suzanne M. Leal, Jing-fei Dong, and Paul F. Bray
Blood, May 2011; 117: 5189 - 5197.

Low-level expression of *let-7a* in gastric cancer and its involvement in tumorigenesis by targeting *RAB40C*

Qiaoyuan Yang, Zhigang Jie, Hong Cao, Anne R. Greenlee, Chengfeng Yang, Fei Zou, and Yiguo Jiang
Carcinogenesis, May 2011; 32: 713 - 722.

IL-23 Receptor Regulation by Let-7f in Human CD4⁺ Memory T Cells

Zhaoxia Li, Feng Wu, Steven R. Brant, and John H. Kwon
J. Immunol., Jun 2011; 186: 6182 - 6190.

Muscle

β -Agonist-associated Reduction in RGS5 Expression Promotes Airway Smooth Muscle Hyper-responsiveness

Zhao Yang, Philip R. Cooper, Gautam Damera, Indranil Mukhopadhyay, Hyeseon Cho, John H. Kehrl, Reynold A. Panettieri, Jr, and Kirk M. Druey
J. Biol. Chem., Apr 2011; 286: 11444 - 11455.

Changes in cross-bridge cycling underlie muscle weakness in patients with tropomyosin 3-based myopathy

Coen A.C. Ottenheijm, Michael W. Lawlor, Ger J.M. Stienen, Henk Granzier, and Alan H. Beggs
Hum. Mol. Genet., May 2011; 20: 2015 - 2025.

Identification of Differentially Regulated Secretome Components During Skeletal Myogenesis

C. Y. X'avia Chan, Olena Masui, Olga Krakovska, Vladimir E. Belozarov, Sebastien Voisin, Shaun Ghanny, Jian Chen, Dharsee Moyez, Peihong Zhu, Kenneth R. Evans, John C. McDermott, and K. W. Michael Siu
Mol. Cell. Proteomics, May 2011; 10: M110.004804.

Myocardin is differentially required for the development of smooth muscle cells and cardiomyocytes

Mark H. Hoofnagle, Ronald L. Nepl, Erica L. Berzin, G. C. Teg Pipes, Eric N. Olson, Brian W. Wamhoff, Avril V. Somlyo, and Gary K. Owens
Am J Physiol Heart Circ Physiol, May 2011; 300: H1707 - H1721.

Tropomyosin variants describe distinct functional subcellular domains in differentiated vascular smooth muscle cells

Cynthia Gallant, Sarah Appel, Philip Graceffa, Paul Leavis, Jim Jung-Ching Lin, Peter W. Gunning, Galina Schevzov, Christine Chaponnier, Jon DeGnore, William Lehman, and Kathleen G. Morgan
Am J Physiol Cell Physiol, Jun 2011; 300: C1356 - C1365.

Rat diaphragm mitochondria have lower intrinsic respiratory rates than mitochondria in limb muscles

Mary L. Garcia-Cazarin, Jorge L. Gamboa, and Francisco H. Andrade
Am J Physiol Regulatory Integrative Comp Physiol, Jun 2011; 300: R1311 - R1315.

Muscle-specific expression of insulin-like growth factor 1 improves outcome in *Lama2^{Dy-w}* mice, a model for congenital muscular dystrophy type 1A

Ajay Kumar, Jenny Yamauchi, Tanya Girgenrath, and Mahasweta Girgenrath
Hum. Mol. Genet., Jun 2011; 20: 2333 - 2343.

Postnatal Changes in the Developing Rat Extraocular Muscles

Carole L. Moncman, Miguel E. Andrade, and Francisco H. Andrade
Invest. Ophthalmol. Vis. Sci., Jun 2011; 52: 3962 - 3969.

Interleukin-13 inhibits proliferation and enhances contractility of human airway smooth muscle cells without change in contractile phenotype

Paul-André Risse, Taisuke Jo, Fernando Suarez, Nobuaki Hirota, Barbara Tolloczko, Pasquale Ferraro, Peter Grutter, and James G. Martin
Am J Physiol Lung Cell Mol Physiol, Jun 2011; 300: L958 - L966.

Neuroscience

The role of Synaptobrevin1/VAMP1 in Ca²⁺-triggered neurotransmitter release at the mouse neuromuscular junction

Yun Liu, Yoshie Sugiura, and Weichun Lin
J. Physiol., Apr 2011; 589: 1603 - 1618.

Altered gene regulation and synaptic morphology in *Drosophila* learning and memory mutants

Zhuo Guan, Lauren K. Buhl, William G. Quinn, and J. Troy Littleton

Learn. Mem., Apr 2011; 18: 191 - 206.**BACE1 Retrograde Trafficking Is Uniquely Regulated by the Cytoplasmic Domain of Sortilin**

Gina M. Finan, Hirokazu Okada, and Tae-Wan Kim

J. Biol. Chem., Apr 2011; 286: 12602 - 12616.**Structural and Functional Analysis of Tomosyn Identifies Domains Important in Exocytotic Regulation**

Antionette L. Williams, Noa Bielopolski, Daphna Meroz, Alice D. Lam, Daniel R. Passmore, Nir Ben-Tal, Stephen A. Ernst, Uri Ashery, and Edward L. Stuenkel

J. Biol. Chem., Apr 2011; 286: 14542 - 14553.**Natural Restoration of Critical Period Plasticity in the Juvenile and Adult Primary Auditory Cortex**

Xiaoming Zhou, Rogerio Panizzutti, Étienne de Villers-Sidani, Caroline Madeira, and Michael M. Merzenich

J. Neurosci., Apr 2011; 31: 5625 - 5634.**Truncation of monocyte chemoattractant protein 1 by plasmin promotes blood–brain barrier disruption**

Yao Yao and Stella E. Tsirka

J. Cell Sci., May 2011; 124: 1486 - 1495.**The Synaptic Proteome during Development and Plasticity of the Mouse Visual Cortex**

Martijn Dahlhaus, Ka Wan Li, Roel C. van der Schors, M. Hadi Saiepour, Pim van Nierop, J. Alexander Heimel,

Josephine M. Hermans, Maarten Loos, August B. Smit, and Christiaan N. Levelt

Mol. Cell. Proteomics, May 2011; 10: M110.005413.**Alzheimer's Disease Brain-Derived Amyloid- β -Mediated Inhibition of LTP *in vivo* Is Prevented by Immunotargeting Cellular Prion Protein**

Andrew E. Barry, Igor Klyubin, Jessica M. Mc Donald, Alexandra J. Mably, Michael A. Farrell, Michael Scott,

Dominic M. Walsh, and Michael J. Rowan

J. Neurosci., May 2011; 31: 7259 - 7263.**Dopamine–Galanin Receptor Heteromers Modulate Cholinergic Neurotransmission in the Rat Ventral Hippocampus**

Estefanía Moreno, Sandra H. Vaz, Ning-Sheng Cai, Carla Ferrada, César Quiroz, Sandeep Kumar Barodia, Nadine Kabbani,

Enric I. Canela, Peter J. McCormick, Carme Lluís, Rafael Franco, Joaquim A. Ribeiro, Ana M. Sebastião, and Sergi Ferré

J. Neurosci., May 2011; 31: 7412 - 7423.**Phosphorylation controls a dual-function polybasic nuclear localization sequence in the adapter protein SH2B1 β to regulate its cellular function and distribution**

Travis J. Maures, Hsiao-Wen Su, Lawrence S. Argetsinger, Sergio Grinstein, and Christin Carter-Su

J. Cell Sci., May 2011; 124: 1542 - 1552.**Long-Term Effects of Repetitive Transcranial Magnetic Stimulation on Markers for Neuroplasticity: Differential Outcomes in Anesthetized and Awake Animals**

Roman Gersner, Elena Kravetz, Jodie Feil, Gaby Pell, and Abraham Zangen

J. Neurosci., May 2011; 31: 7521 - 7526.**Soluble A β Oligomers Inhibit Long-Term Potentiation through a Mechanism Involving Excessive Activation of Extrasynaptic NR2B-Containing NMDA Receptors**

Shaomin Li, Ming Jin, Thomas Koeglsperger, Nina E. Shepardson, Ganesh M. Shankar, and Dennis J. Selkoe

J. Neurosci., May 2011; 31: 6627 - 6638.**Overexpression of the Dominant-Negative Form of Interferon Regulatory Factor 1 in Oligodendrocytes Protects against Experimental Autoimmune Encephalomyelitis**

Zihua Ren, Yan Wang, Duan Tao, David Liebenson, Thomas Liggett, Rajendra Goswami, Robert Clarke, Dusan Stefanoski,

and Roumen Balabanov

J. Neurosci., Jun 2011; 31: 8329 - 8341.

ApoE isoform-specific regulation of regeneration in the peripheral nervous system

Laura H. Comley, Heidi R. Fuller, Thomas M. Wishart, Chantal A. Mutsaers, Derek Thomson, Ann K. Wright, Richard R. Ribchester, Glenn E. Morris, Simon H. Parson, Karen Horsburgh, and Thomas H. Gillingwater
Hum. Mol. Genet., Jun 2011; 20: 2406 - 2421.

Scapinin-induced Inhibition of Axon Elongation Is Attenuated by Phosphorylation and Translocation to the Cytoplasm

Hovik Farghaian, Yu Chen, Ada W. Y. Fu, Amy K. Y. Fu, Jacques P. K. Ip, Nancy Y. Ip, Ann M. Turnley, and Adam R. Cole
J. Biol. Chem., Jun 2011; 286: 19724 - 19734.

Specific Regulation of NRG1 Isoform Expression by Neuronal Activity

Xihui Liu, Ryan Bates, Dong-Min Yin, Chengyong Shen, Fay Wang, Nan Su, Sergei A. Kirov, Yuling Luo, Jian-Zhi Wang, Wen-Cheng Xiong, and Lin Mei
J. Neurosci., Jun 2011; 31: 8491 - 8501.

Wild-type HTT modulates the enzymatic activity of the neuronal palmitoyl transferase HIP14

Kun Huang, Shaun S. Sanders, Rujun Kang, Jeffrey B. Carroll, Liza Sutton, Junmei Wan, Roshni Singaraja, Fiona B. Young, Lili Liu, Alaa El-Husseini, Nicholas G. Davis, and Michael R. Hayden
Hum. Mol. Genet., Jun 2011; 10.1093/hmg/ddr242.

Nephrology/Kidney

TGF- β 1-activated kinase-1 regulates inflammation and fibrosis in the obstructed kidney

Frank Y. Ma, Greg H. Tesch, Elyce Ozols, Min Xie, Michael D. Schneider, and David J. Nikolic-Paterson
Am J Physiol Renal Physiol, Jun 2011; 300: F1410 - F1421.

Phosphatases

Sts-2 Is a Phosphatase That Negatively Regulates Zeta-associated Protein (ZAP)-70 and T Cell Receptor Signaling Pathways

Boris San Luis, Ben Sondgeroth, Nicolas Nassar, and Nick Carpino
J. Biol. Chem., May 2011; 286: 15943 - 15954.

The E3 Ubiquitin Ligase- and Protein Phosphatase 2A (PP2A)-binding Domains of the Alpha4 Protein Are Both Required for Alpha4 to Inhibit PP2A Degradation

Michele LeNoue-Newton, Guy R. Watkins, Ping Zou, Katherine L. Germane, Lisa R. McCorvey, Brian E. Wadzinski, and Benjamin W. Spiller
J. Biol. Chem., May 2011; 286: 17665 - 17671.

Phosphatase-Resistant Gap Junctions Inhibit Pathological Remodeling and Prevent Arrhythmias

Benjamin F. Remo, Jiaxiang Qu, Frank M. Volpicelli, Steven Giovannone, Daniel Shin, Joshua Lader, Fang-yu Liu, Jie Zhang, Danielle S. Lent, Gregory E. Morley, and Glenn I. Fishman
Circ. Res., Jun 2011; 108: 1459 - 1466.

LipA, a Tyrosine and Lipid Phosphatase Involved in the Virulence of *Listeria monocytogenes*

Renate Kastner, Olivier Dussurget, Cristel Archambaud, Elisabeth Kernbauer, Didier Soulat, Pascale Cossart, and Thomas Decker
Infect. Immun., Jun 2011; 79: 2489 - 2498.

Regulation of Src Family Kinases Involved in T Cell Receptor Signaling by Protein-tyrosine Phosphatase CD148

Ondrej Stepanek, Tomas Kalina, Peter Draber, Tereza Skopцова, Karel Svojgr, Pavla Angelisova, Vaclav Horejsi, Arthur Weiss, and Tomas Brdicka
J. Biol. Chem., Jun 2011; 286: 22101 - 22112.

Phosphatase-Resistant Gap Junctions Inhibit Pathological Remodeling and Prevent Arrhythmias

Benjamin F. Remo, Jiaxiang Qu, Frank M. Volpicelli, Steven Giovannone, Daniel Shin, Joshua Lader, Fang-yu Liu, Jie Zhang, Danielle S. Lent, Gregory E. Morley, and Glenn I. Fishman
Circ. Res., Jun 2011; 108: 1459 - 1466.

Receptors

Targeted Therapy of VEGFR2 and EGFR Significantly Inhibits Growth of Anaplastic Thyroid Cancer in an Orthotopic Murine Model

Maria K. Gule, Yunyun Chen, Daisuke Sano, Mitchell J. Frederick, Ge Zhou, Mei Zhao, Zvonimir L. Milas, Chad E. Galer, Ying C. Henderson, Samar A. Jasser, David L. Schwartz, James A. Bankson, Jeffrey N. Myers, and Stephen Y. Lai
Clin. Cancer Res., Apr 2011; 17: 2281 - 2291.

Differential responses of epithelial Sertoli cells of the rat testis to Toll-like receptor 2 and 4 ligands: Implications for studies of testicular inflammation using bacterial lipopolysaccharides

Wendy R. Winnall, Julie A. Muir, and Mark P. Hedger
Innate Immunity, Apr 2011; 17: 123 - 136.

KISS1R Intracellular Trafficking and Degradation: Effect of the Arg386Pro Disease-Associated Mutation

Suzy D. C. Bianco, Lauren Vandepas, Mayrin Correa-Medina, Balázs Gereben, Abir Mukherjee, Wendy Kuohung, Rona Carroll, Milena G. Teles, Ana Claudia Latronico, and Ursula B. Kaiser
Endocrinology, Apr 2011; 152: 1616 - 1626.

Soluble Form of the (Pro)Renin Receptor Is Augmented in the Collecting Duct and Urine of Chronic Angiotensin II-Dependent Hypertensive Rats

Alexis A. Gonzalez, Lucienne S. Lara, Christina Luffman, Dale M. Seth, and Minolfa C. Prieto
Hypertension, Apr 2011; 57: 859 - 864.

Agonist-Biased Signaling at the Histamine H₄ Receptor: JNJ7777120 Recruits β -Arrestin without Activating G Proteins

Elizabeth M. Rosethorne and Steven J. Charlton
Mol. Pharmacol., Apr 2011; 79: 749 - 757.

ARH cooperates with AP-1B in the exocytosis of LDLR in polarized epithelial cells

Richard S. Kang and Heike Fölsch
J. Cell Biol., Apr 2011; 193: 51 - 60.

Urocortin 1 modulates immunosignaling in a rat model of colitis via corticotropin-releasing factor receptor 2

Jen Chang, Melanie R. Adams, Matthew S. Clifton, Min Liao, Julia H. Brooks, Burcu Hasdemir, and Aditi Bhargava
Am J Physiol Gastrointest Liver Physiol, May 2011; 300: G884 - G894.

Calcium/Calmodulin-Dependent Protein Kinase II Mediates Group I Metabotropic Glutamate Receptor-Dependent Protein Synthesis and Long-Term Depression in Rat Hippocampus

Bruce G. Mockett, Diane Guévremont, Magdalena Wutte, Sarah R. Hulme, Joanna M. Williams, and Wickliffe C. Abraham
J. Neurosci., May 2011; 31: 7380 - 7391.

A Multi-domain Fragment of Nogo-A Protein Is a Potent Inhibitor of Cortical Axon Regeneration via Nogo Receptor 1

Eric A. Huebner, Byung G. Kim, Philip J. Duffy, Rebecca H. Brown, and Stephen M. Strittmatter
J. Biol. Chem., May 2011; 286: 18026 - 18036.

IL-23 Receptor Regulation by Let-7f in Human CD4⁺ Memory T Cells

Zhaoxia Li, Feng Wu, Steven R. Brant, and John H. Kwon
J. Immunol., Jun 2011; 186: 6182 - 6190.

Antitumor and Antiangiogenic Activities of BMS-690514, an Inhibitor of Human EGF and VEGF Receptor Kinase Families

Tai W. Wong, Francis Y. Lee, Stuart Emanuel, Craig Fairchild, Joseph Fargnoli, Brian Fink, Ashvinikumar Gavai, Amy Hammell, Benjamin Henley, Christine Hilt, John T. Hunt, Bala Krishnan, Daniel Kukral, Anne Lewin, Harold Malone, Derek Norris, Simone Oppenheimer, Gregory Vite, and Chiang Yu
Clin. Cancer Res., Jun 2011; 17: 4031 - 4041.

Pharmacological inhibition of sphingosine kinase isoforms alters estrogen receptor signaling in human breast cancer

James W Antoon, William D Meacham, Melyssa R Bratton, Evelyn M Slaughter, Lyndsay V Rhodes, Hasina B Ashe, Thomas E Wiese, Matthew E Burow, and Barbara S Beckman
J. Mol. Endocrinol., Jun 2011; 46: 205 - 216.

Kinase Suppressor of Ras 1 (KSR1) Regulates PGC1 α and Estrogen-Related Receptor α To Promote Oncogenic Ras-Dependent Anchorage-Independent Growth

Kurt W. Fisher, Binita Das, Robert L. Kortum, Oleg V. Chaika, and Robert E. Lewis
Mol. Cell. Biol., Jun 2011; 31: 2453 - 2461.

Regulation of Src Family Kinases Involved in T Cell Receptor Signaling by Protein-tyrosine Phosphatase CD148

Ondrej Stepanek, Tomas Kalina, Peter Draber, Tereza Skopcova, Karel Svojgr, Pavla Angelisova, Vaclav Horejsi, Arthur Weiss, and Tomas Brdicka
J. Biol. Chem., Jun 2011; 286: 22101 - 22112.

RNAi

MicroRNA-29c Is a Signature MicroRNA under High Glucose Conditions That Targets Sprouty Homolog 1, and its *in vivo* Knockdown Prevents Progression of Diabetic Nephropathy

Jianyin Long, Yin Wang, Wenjian Wang, Benny H. J. Chang, and Farhad R. Danesh
J. Biol. Chem., Apr 2011; 286: 11837 - 11848.

Depletion of PINK1 affects mitochondrial metabolism, calcium homeostasis and energy maintenance

Bavo Heeman, Chris Van den Haute, Sarah-Ann Aelvoet, Federica Valsecchi, Richard J. Rodenburg, Veerle Reumers, Zeger Debyser, Geert Callewaert, Werner J. H. Koopman, Peter H. G. M. Willems, and Veerle Baekelandt
J. Cell Sci., Apr 2011; 124: 1115 - 1125.

BACE1 Retrograde Trafficking Is Uniquely Regulated by the Cytoplasmic Domain of Sortilin

Gina M. Finan, Hirokazu Okada, and Tae-Wan Kim
J. Biol. Chem., Apr 2011; 286: 12602 - 12616.

Platelet Protein Kinase C- ϕ Deficiency With Human RUNX1 Mutation: *PRKCQ* Is a Transcriptional Target of RUNX1

Gauthami Jalagadugula, Guangfen Mao, Gurpreet Kaur, Danny N. Dhanasekaran, and A. Koneti Rao
Arterioscler Thromb Vasc Biol, Apr 2011; 31: 921 - 927.

Notch1-mediated Signaling Induces MHC Class II Expression through Activation of Class II Transactivator Promoter III in Mast Cells

Nobuhiro Nakano, Chiharu Nishiyama, Hideo Yagita, Akemi Koyanagi, Hideoki Ogawa, and Ko Okumura
J. Biol. Chem., Apr 2011; 286: 12042 - 12048.

Role of Inflammation and Insulin Resistance in Endothelial Progenitor Cell Dysfunction

Cyrus V. Desouza, Frederick G. Hamel, Keshore Bidasee, and Kelly O'Connell
Diabetes, Apr 2011; 60: 1286 - 1294.

Inhibition of renal caveolin-1 reduces natriuresis and produces hypertension in sodium-loaded rats

John J. Gildea, Brandon A. Kemp, Nancy L. Howell, Robert E. Van Sciver, Robert M. Carey, and Robin A. Felder
Am J Physiol Renal Physiol, Apr 2011; 300: F914 - F920.

Antisense oligonucleotide reduction of apoB-ameliorated atherosclerosis in LDL receptor-deficient mice

Adam E. Mullick, Wuxia Fu, Mark J. Graham, Richard G. Lee, Donna Witchell, Thomas A. Bell, Charles P. Whipple, and Rosanne M. Crooke
J. Lipid Res., May 2011; 52: 885 - 896.

Tumour suppressor ING1b maintains genomic stability upon replication stress

Ronald P. C. Wong, Hanyang Lin, Shahram Khosravi, Brad Piche, Seyed Mehdi Jafarnejad, David W. C. Chen, and Gang Li
Nucleic Acids Res., May 2011; 39: 3632 - 3642.

Mechanisms of single-stranded phosphorothioate modified antisense oligonucleotide accumulation in hepatocytes

Erich Koller, Thomas M. Vincent, Alfred Chappell, Soma De, Muthiah Manoharan, and C. Frank Bennett
Nucleic Acids Res., Jun 2011; 39: 4795 - 4807.

Intracellular targeting of peroxiredoxin 6 to lysosomal organelles requires MAPK activity and binding to 14-3-3ε

Elena M. Sorokina, Sheldon I. Feinstein, Suiping Zhou, and Aron B. Fisher
Am J Physiol Cell Physiol, Jun 2011; 300: C1430 - C1441.

Wild-type HTT modulates the enzymatic activity of the neuronal palmitoyl transferase HIP14

Kun Huang, Shaun S. Sanders, Rujun Kang, Jeffrey B. Carroll, Liza Sutton, Junmei Wan, Roshni Singaraja, Fiona B. Young, Lili Liu, Alaa El-Husseini, Nicholas G. Davis, and Michael R. Hayden
Hum. Mol. Genet., Jun 2011; 10.1093/hmg/ddr242.

HOXC9 Links Cell-Cycle Exit and Neuronal Differentiation and Is a Prognostic Marker in Neuroblastoma

Ling Mao, Jane Ding, Yunhong Zha, Liqun Yang, Brian A. McCarthy, William King, Hongjuan Cui, and Han-Fei Ding
Cancer Res., Jun 2011; 71: 4314 - 4324.

Expanding the Paradigm for Estrogen Receptor Binding and Transcriptional Activation

S. R. Joshi, R. B. Ghattamaneni, and W. M. Scovell
Mol. Endocrinol., Jun 2011; 25: 980 - 994.

Distinct roles of cadherin-6 and E-cadherin in tubulogenesis and lumen formation

Liwei Jia, Fengming Liu, Steen H. Hansen, Martin B.A. ter Beest, and Mirjam M.P. Zegers
Mol. Biol. Cell, Jun 2011; 22: 2031 - 2041.

Transcription Factors

Notch1-mediated Signaling Induces MHC Class II Expression through Activation of Class II Transactivator Promoter III in Mast Cells

Nobuhiro Nakano, Chiharu Nishiyama, Hideo Yagita, Akemi Koyanagi, Hideoki Ogawa, and Ko Okumura
J. Biol. Chem., Apr 2011; 286: 12042 - 12048.

C/EBPβ Mediates Growth Hormone-Regulated Expression of Multiple Target Genes

Tracy X. Cui, Grace Lin, Christopher R. LaPensee, Anda-Alexandra Calinescu, Maanjot Rathore, Cale Streeter, Graciela Piwien-Pilipuk, Nathan Lanning, Hui Jin, Christin Carter-Su, Zhaohui S. Qin, and Jessica Schwartz
Mol. Endocrinol., Apr 2011; 25: 681 - 693.

STAT3 Mediates Resistance to MEK Inhibitor through MicroRNA miR-17

Bingbing Dai, Jieru Meng, Michael Peyton, Luc Girard, William G. Bornmann, Lin Ji, John D. Minna, Bingliang Fang, and Jack A. Roth
Cancer Res., May 2011; 71: 3658 - 3668.

Stat3 signaling in acute myeloid leukemia: ligand-dependent and -independent activation and induction of apoptosis by a novel small-molecule Stat3 inhibitor

Michele S. Redell, Marcos J. Ruiz, Todd A. Alonzo, Robert B. Gerbing, and David J. Tweardy
Blood, May 2011; 117: 5701 - 5709.

Smad3 Dosage Determines Androgen Responsiveness and Sets the Pace of Postnatal Testis Development

Catherine Itman, Chin Wong, Briony Hunyadi, Matthias Ernst, David A. Jans, and Kate L. Loveland
Endocrinology, May 2011; 152: 2076 - 2089.

Cytokine-induced Paracrystals Prolong the Activity of Signal Transducers and Activators of Transcription (STAT) and Provide a Model for the Regulation of Protein Solubility by Small Ubiquitin-like Modifier (SUMO)

Mathias Droscher, Andreas Begitt, Andreas Marg, Martin Zacharias, and Uwe Vinkemeier
J. Biol. Chem., May 2011; 286: 18731 - 18746.

Suberoylanilide Hydroxamic Acid (Vorinostat) Up-regulates Progranulin Transcription:**RATIONAL THERAPEUTIC APPROACH TO FRONTOTEMPORAL DEMENTIA**

Basar Cenik, Chantelle F. Sephton, Colleen M. Dewey, Xunde Xian, Shuguang Wei, Kimberley Yu, Wenze Niu, Giovanni Coppola, Sarah E. Coughlin, Suzee E. Lee, Daniel R. Dries, Sandra Almeida, Daniel H. Geschwind, Fen-Biao Gao, Bruce L. Miller, Robert V. Farese, Jr, Bruce A. Posner, Gang Yu, and Joachim Herz
J. Biol. Chem., May 2011; 286: 16101 - 16108.

AMP-activated Protein Kinase Regulates β -Catenin Transcription via Histone Deacetylase 5

Jun-Xing Zhao, Wan-Fu Yue, Mei-Jun Zhu, and Min Du
J. Biol. Chem., May 2011; 286: 16426 - 16434.

Mapping the κ B Kinase β (IKK β)-binding Interface of the B14 Protein, a Vaccinia Virus Inhibitor of IKK β -mediated Activation of Nuclear Factor κ B

Camilla T. O. Benfield, Daniel S. Mansur, Laura E. McCoy, Brian J. Ferguson, Mohammad W. Bahar, Asa P. Oldring, Jonathan M. Grimes, David I. Stuart, Stephen C. Graham, and Geoffrey L. Smith
J. Biol. Chem., Jun 2011; 286: 20727 - 20735.

Epithelial Cell Organization Suppresses Myc Function by Attenuating Myc Expression

David R. Simpson, Min Yu, Siyuan Zheng, Zhongming Zhao, Senthil K. Muthuswamy, and William P. Tansey
Cancer Res., Jun 2011; 71: 3822 - 3830.

Quantification of Activated NF- κ B/RelA Complexes Using ssDNA Aptamer Affinity – Stable Isotope Dilution – Selected Reaction Monitoring – Mass Spectrometry

Yingxin Zhao, Steven G. Widen, Mohammad Jamaluddin, Bing Tian, Thomas G. Wood, Chukwudi B. Edeh, and Allan R. Brasier
Mol. Cell. Proteomics, Jun 2011; 10: M111.008771.

The Pleckstrin Homology Domain in the SKAP55 Adapter Protein Defines the Ability of the Adapter Protein ADAP To Regulate Integrin Function and NF- κ B Activation

Brandon J. Burbach, Rupa Srivastava, Melissa A. Ingram, Jason S. Mitchell, and Yoji Shimizu
J. Immunol., Jun 2011; 186: 6227 - 6237.

p53 Negatively Regulates RGS13 Protein Expression in Immune Cells

Shoko Iwaki, Yunbiao Lu, Zhihui Xie, and Kirk M. Druey
J. Biol. Chem., Jun 2011; 286: 22219 - 22226.

Cardiac-specific mindin overexpression attenuates cardiac hypertrophy via blocking AKT/GSK3 β and TGF- β 1-Smad signalling

Ling Yan, Xiang Wei, Qi-Zhu Tang, Jinghua Feng, Yan Zhang, Chen Liu, Zhou-Yan Bian, Lian-Feng Zhang, Manyin Chen, Xue Bai, Ai-Bing Wang, John Fassett, Yingjie Chen, You-Wen He, Qinglin Yang, Peter P. Liu, and Hongliang Li
Cardiovasc Res, Jun 2011; 10.1093/cvr/cvr159.

TGF- β 1-activated kinase-1 regulates inflammation and fibrosis in the obstructed kidney

Frank Y. Ma, Greg H. Tesch, Elyce Ozols, Min Xie, Michael D. Schneider, and David J. Nikolic-Paterson
Am J Physiol Renal Physiol, Jun 2011; 300: F1410 - F1421.

The *Agrobacterium tumefaciens* Transcription Factor B1cR Is Regulated via Oligomerization

Yi Pan, Valena Fiscus, Wuyi Meng, Zhida Zheng, Lian-Hui Zhang, Clay Fuqua, and Lingling Chen
J. Biol. Chem., Jun 2011; 286: 20431 - 20440.

Tight Control of STAT5 Activity Determines Human CD34-Derived Interstitial Dendritic Cell and Langerhans Cell Development

Lianne van de Laar, Aniek van den Bosch, Albertus T. J. Wierenga, Harry L. A. Janssen, Paul J. Coffey, and Andrea M. Woltman
J. Immunol., Jun 2011; 186: 7016 - 7024.

Mutations in the *NKX2.5* Gene and the *PAX8* Promoter in a Girl with Thyroid Dysgenesis

Pia Hermanns, Helmut Grasberger, Samuel Refetoff, and Joachim Pohlenz
J. Clin. Endocrinol. Metab., Jun 2011; 96: E977 - E981.

Ubiquitination and Sumoylation

Structural and Functional Analysis of Tomosyn Identifies Domains Important in Exocytotic Regulation

Antionette L. Williams, Noa Bielopolski, Daphna Meroz, Alice D. Lam, Daniel R. Passmore, Nir Ben-Tal, Stephen A. Ernst, Uri Ashery, and Edward L. Stuenkel
J. Biol. Chem., Apr 2011; 286: 14542 - 14553.

The E3 Ubiquitin Ligase- and Protein Phosphatase 2A (PP2A)-binding Domains of the Alpha4 Protein Are Both Required for Alpha4 to Inhibit PP2A Degradation

Michele LeNoue-Newton, Guy R. Watkins, Ping Zou, Katherine L. Germane, Lisa R. McCorvey, Brian E. Wadzinski, and Benjamin W. Spiller
J. Biol. Chem., May 2011; 286: 17665 - 17671.

Cytokine-induced Paracrystals Prolong the Activity of Signal Transducers and Activators of Transcription (STAT) and Provide a Model for the Regulation of Protein Solubility by Small Ubiquitin-like Modifier (SUMO)

Mathias Droscher, Andreas Begitt, Andreas Marg, Martin Zacharias, and Uwe Vinkemeier
J. Biol. Chem., May 2011; 286: 18731 - 18746.

Activities of ICP0 Involved in the Reversal of Silencing of Quiescent Herpes Simplex Virus 1

Michael W. Ferenczy, Daniel J. Ranayhossaini, and Neal A. DeLuca
J. Virol., May 2011; 85: 4993 - 5002.

Sequential requirements for the GTPase domain of the mitofusin Fzo1 and the ubiquitin ligase SCF^{Mdm30} in mitochondrial outer membrane fusion

Mickael M. Cohen, Elizabeth A. Amiot, Adam R. Day, Guillaume P. Leboucher, Erin N. Pryce, Michael H. Glickman, J. Michael McCaffery, Janet M. Shaw, and Allan M. Weissman
J. Cell Sci., May 2011; 124: 1403 - 1410.

Impairment of Human Immunodeficiency Virus Type-1 Integrase SUMOylation Correlates with an Early Replication Defect

Alessia Zamborlini, Audrey Coiffic, Guillaume Beauclair, Olivier Delelis, Joris Paris, Yashuiro Koh, Fabian Magne, Marie-Lou Giron, Joelle Tobaly-Tapiero, Eric Deprez, Stephane Emiliani, Alan Engelman, Hugues deThé, and Ali Saïb
J. Biol. Chem., Jun 2011; 286: 21013 - 21022.

Prefoldin Subunits Are Protected from Ubiquitin-Proteasome System-mediated Degradation by Forming Complex with Other Constituent Subunits

Makoto Miyazawa, Erika Tashiro, Hirotake Kitaura, Hiroshi Maita, Hiroo Suto, Sanae M. M. Iguchi-Ariga, and Hiroyoshi Ariga
J. Biol. Chem., Jun 2011; 286: 19191 - 19203.

Virology/Microbiology

Structural basis for the neutralization and genotype specificity of hepatitis E virus

Xuhua Tang, Chunyan Yang, Ying Gu, Cuiling Song, Xiao Zhang, Yingbin Wang, Jun Zhang, Choy Leong Hew, Shaowei Li, Ningshao Xia, and J. Sivaraman
PNAS, Jun 2011; 108: 10266 - 10271.

Multiple Interactions between the ESCRT Machinery and Arrestin-Related Proteins: Implications for PPXY-Dependent Budding

Susanne Rauch and Juan Martin-Serrano
J. Virol., Apr 2011; 85: 3546 - 3556.

Hepatitis B Virus X Protein Enhances Cisplatin-Induced Hepatotoxicity via a Mechanism Involving Degradation of Mcl-1

Liang Hu, Lei Chen, Liang Li, HanYong Sun, GuangZhen Yang, YanXin Chang, QianQian Tu, MengChao Wu, and HongYang Wang
J. Virol., Apr 2011; 85: 3214 - 3228.

IL-15 and Type I Interferon Are Required for Activation of Tumoricidal NK Cells by Virus-Infected Dendritic Cells

Jeanette E. Boudreau, Kyle B. Stephenson, Fuan Wang, Ali A. Ashkar, Karen L. Mossman, Laurel L. Lenz, Kenneth L. Rosenthal, Jonathan L. Bramson, Brian D. Lichty, and Yonghong Wan
Cancer Res., Apr 2011; 71: 2497 - 2506.

LIM Kinase 1 Modulates Cortical Actin and CXCR4 Cycling and Is Activated by HIV-1 to Initiate Viral Infection

Paul J. Vorster, Jia Guo, Alyson Yoder, Weifeng Wang, Yanfang Zheng, Xuehua Xu, Dongyang Yu, Mark Spear, and Yuntao Wu
J. Biol. Chem., Apr 2011; 286: 12554 - 12564.

C/EBP β Mediates Growth Hormone-Regulated Expression of Multiple Target Genes

Tracy X. Cui, Grace Lin, Christopher R. LaPensee, Anda-Alexandra Calinescu, Maanjot Rathore, Cale Streeter, Graciela Pivien-Pilipuk, Nathan Lanning, Hui Jin, Christin Carter-Su, Zhaohui S. Qin, and Jessica Schwartz
Mol. Endocrinol., Apr 2011; 25: 681 - 693.

Activities of ICP0 Involved in the Reversal of Silencing of Quiescent Herpes Simplex Virus 1

Michael W. Ferenczy, Daniel J. Ranayhossaini, and Neal A. DeLuca
J. Virol., May 2011; 85: 4993 - 5002.

Inactivation of hantaviruses by *N*-ethylmaleimide preserves virion integrity

Tomas Strandin, Jussi Hepojoki, Hao Wang, Antti Vaeheri, and Hilikka Lankinen
J. Gen. Virol., May 2011; 92: 1189 - 1198.

Residues of the UL25 Protein of Herpes Simplex Virus That Are Required for Its Stable Interaction with Capsids

Shelley K. Cockrell, Jamie B. Huffman, Katerina Toropova, James F. Conway, and Fred L. Homa
J. Virol., May 2011; 85: 4875 - 4887.

A Rab11- and Microtubule-Dependent Mechanism for Cytoplasmic Transport of Influenza A Virus Viral RNA

Maria Joao Amorim, Emily A. Bruce, Eliot K. C. Read, Ágnes Foeglein, Robert Mahen, Amanda D. Stuart, and Paul Digard
J. Virol., May 2011; 85: 4143 - 4156.

Formation of a Stable RuvA Protein Double Tetramer Is Required for Efficient Branch Migration *in vitro* and for Replication Fork Reversal *in vivo*

Alison S. Bradley, Zeynep Baharoglu, Andrew Niewiarowski, Bénédicte Michel, and Irina R. Tsaneva
J. Biol. Chem., Jun 2011; 286: 22372 - 22383.

DNA Repair and Global Sumoylation Are Regulated by Distinct Ubc9 Noncovalent Complexes

John Prudden, J. Jefferson P. Perry, Minghua Nie, Ajay A. Vashisht, Andrew S. Arvai, Chiharu Hitomi, Grant Guenther, James A. Wohlschlegel, John A. Tainer, and Michael N. Boddy
Mol. Cell. Biol., Jun 2011; 31: 2299 - 2310.

The Herpes Simplex Virus 1 vhs Protein Enhances Translation of Viral True Late mRNAs and Virus Production in a Cell Type-Dependent Manner

Bianca Dauber, Jerry Pelletier, and James R. Smiley
J. Virol., Jun 2011; 85: 5363 - 5373.

Overlapping signals for translational regulation and packaging of influenza A virus segment 2

Helen M. Wise, Cyril Barbezange, Brett W. Jagger, Rosa M. Dalton, Julia R. Gog, Martin D. Curran, Jeffery K. Taubenberger, Emma C. Anderson, and Paul Digard
Nucleic Acids Res., Jun 2011; 10.1093/nar/gkr487.

Rescue of Infectious Particles from Preassembled Alphavirus Nucleocapsid Cores

Jonathan E. Snyder, Odisse Azizgolshani, Bingbing Wu, Yingpei He, Aih Cheun Lee, Joyce Jose, Daniel M. Suter, Charles M. Knobler, William M. Gelbart, and Richard J. Kuhn
J. Virol., Jun 2011; 85: 5773 - 5781.

Characterization of Empty Adenovirus Particles Assembled in the Absence of a Functional Adenovirus IVa2 Protein

Philomena Ostapchuk, Matthew Almond, and Patrick Hearing
J. Virol., Jun 2011; 85: 5524 - 5531.

APPLICATIONS

In-Cell™ Western Assays

Role of Inflammation and Insulin Resistance in Endothelial Progenitor Cell Dysfunction

Cyrus V. Desouza, Frederick G. Hamel, Keshore Bidasee, and Kelly O'Connell
Diabetes, Apr 2011; 60: 1286 - 1294.

The Protease-activated Receptor-2-specific Agonists 2-Aminothiazol-4-yl-LIGRL-NH₂ and 6-Aminonicotinyl-LIGRL-NH₂ Stimulate Multiple Signaling Pathways to Induce Physiological Responses *in vitro* and *in vivo*

Andrea N. Flynn, Dipti V. Tillu, Marina N. Asiedu, Justin Hoffman, Josef Vagner, Theodore J. Price, and Scott Boitano
J. Biol. Chem., May 2011; 286: 19076 - 19088.

Multipathway Model Enables Prediction of Kinase Inhibitor Cross-Talk Effects on Migration of Her2-Overexpressing Mammary Epithelial Cells

Kumar N, Afeyan R, Kim HD, Lauffenburger DA.
Molecular Pharmacology 73:1668–1678 (2008).

IRDye® 800CW Toxicity Study

Single-Dose Intravenous Toxicity Study of IRDye 800CW in Sprague-Dawley Rats.

MV Marshall, D Draney, EM Sevick-Muraca, and DM Olive
Mol Imaging Biol, Apr 2010.

Membrane Stripping

Optimization of stripping/reprobing for LI-COR® Western blot analysis

Michael C McCauley, Ari Nalbandian, Gabby Brum, Shan Lu, and Yinsheng Wan
FASEB J, Apr 2010; 24: 703.2.

Micro-Western Assay (alternative to reverse phase arrays)

Systems analysis of EGF receptor signaling dynamics with microwestern arrays.

MF Ciaccio, JP Wagner, CP Chuu, DA Lauffenburger, and RB Jones
Nat Methods, Feb 2010; 7(2): 148-55.

MousePod® Accessories

High Plasma Levels and Effective Lymphatic Uptake of Docetaxel in an Orally Available Nanotransporter Formulation

Taher Nassar, Suha Attili-Qadri, Oshrat Harush-Frenkel, Shimon Farber, Shimon Lecht, Philip Lazarovici, and Simon Benita
Cancer Res., Apr 2011; 71: 3018 - 3028.

Functional divergence between 2 chemokines is conferred by single amino acid change

Alexandre Dubrac, Cathy Quemener, Eric Lacazette, Frédéric Lopez, Catherine Zanibellato, Wen-Guey Wu, Andr as Bikfalvi, and Herv  Prats
Blood, Nov 2010; 116: 4703 - 4711.

Antitumor Activity of Gemcitabine and Oxaliplatin Is Augmented by Thymoquinone in Pancreatic Cancer

Sanjeev Banerjee, Ahmed O. Kaseb, Zhiwei Wang, Deujan Kong, Mussop Mohammad, Subhash Padhye, Fazlul H. Sarkar, and Ramzi M. Mohammad
Cancer Res 69: 5575 (2009)

Reverse Phase Protein Arrays

Lysate Microarrays Enable High-throughput, Quantitative Investigations of Cellular Signaling

Mark Sevecka, Alejandro Wolf-Yadlin, and Gavin MacBeath

Mol. Cell. Proteomics, Apr 2011; 10: M110.005363.

Systems Biology Reveals New Strategies for Personalizing Cancer Medicine and Confirms the Role of PTEN in Resistance to Trastuzumab

Dana Faratian, Alexey Goltsov, Galina Lebedeva, Anatoly Sorokin, Stuart Moodie, Peter Mullen, Charlene Kay, In Hwa Um, Simon Langdon, Igor Goryanin, and David J. Harrison

Cancer Res., Aug 2009; 69: 6713 - 6720.

RNA/Protein Interactions

Near-infrared (NIR) dye-labeled RNAs identify binding of ZBP1 to the noncoding Y3-RNA

Marcel Köhn, Marcell Lederer, Kristin Wächter, and Stefan Hüttelmaier

RNA, Jul 2010; 16: 1420 - 1428.

siRNA Screening

Quantitative evaluation of siRNA delivery *in vivo*

Yi Pei, Paula J. Hancock, Hangchun Zhang, René Bartz, Craig Cherrin, Nathalie Innocent, Colin J. Pomerantz, Jessica Seitzer, Martin L. Koser, Marc T. Abrams, Yan Xu, Nelly A. Kuklin, Paul A. Burke, Alan B. Sachs, Laura Sepp-Lorenzino, and Stanley F. Barnett

RNA, Dec 2010; 16: 2553 - 2563.

Western (Quantitative, Molecules/Cell)

Quantitative analyses reveal the importance of regulated Hdmx degradation for P53 activation

Yunyuan V. Wang, Mark Wade, EeTsin Wong, Yao-Cheng Li, Luo Wei Rodewald, and Geoffrey M. Wahl

PNAS, Jul 2007; 104: 12365 – 12370.

PNAS and Science Publications for the quarter (7)

GENETICS

Acidic nuclear phosphoprotein 32kDa (ANP32) β -deficient mouse reveals a hierarchy of ANP32 importance in mammalian development

Patrick T. Reilly, Samia Afzal, Chiara Gorrini, Koren Lui, Yury V. Bukhman, Andrew Wakeham, Jillian Haight, Teo Wei Ling, Carol C. Cheung, Andrew J. Elia, Patricia V. Turner, and Tak Wah Mak
PNAS, Jun 2011; 108: 10243 - 10248.

MEDICAL SCIENCES

Mutations in lipoprotein lipase that block binding to the endothelial cell transporter GPIHBP1

Constance V. Voss, Brandon S. J. Davies, Shelly Tat, Peter Gin, Loren G. Fong, Christopher Pelletier, Charlene D. Mottler, André Bensadoun, Anne P. Beigneux, and Stephen G. Young
PNAS, May 2011; 108: 7980 - 7984.

***In vivo* diagnosis of murine pancreatic intraepithelial neoplasia and early-stage pancreatic cancer by molecular imaging**

Stefan Eser, Marlena Messer, Philipp Eser, Alexander von Werder, Barbara Seidler, Monther Bajbouj, Roger Vogelmann, Alexander Meining, Johannes von Burstin, Hana Algül, Philipp Pagel, Angelika E. Schnieke, Irene Esposito, Roland M. Schmid, Günter Schneider, and Dieter Saur
PNAS, Jun 2011; 108: 9945 - 9950.

Structural basis for the neutralization and genotype specificity of hepatitis E virus

Xuhua Tang, Chunyan Yang, Ying Gu, Cuiling Song, Xiao Zhang, Yingbin Wang, Jun Zhang, Choy Leong Hew, Shaowei Li, Ningshao Xia, and J. Sivaraman
PNAS, Jun 2011; 108: 10266 - 10271.

NEUROSCIENCE

Impaired NMDA receptor transmission alters striatal synapses and DISC1 protein in an age-dependent manner

Amy J. Ramsey, Marija Milenkovic, Ana F. Oliveira, Yasmin Escobedo-Lozoya, Saurav Seshadri, Ali Salahpour, Akira Sawa, Ryohei Yasuda, and Marc G. Caron
PNAS, Apr 2011; 108: 5795 - 5800.

RAS-converting enzyme 1-mediated endoproteolysis is required for trafficking of rod phosphodiesterase 6 to photoreceptor outer segments

Jeffrey R. Christiansen, Saravanan Koldaivelu, Martin O. Bergo, and Visvanathan Ramamurthy
PNAS, May 2011; 108: 8862 - 8866.

PHYSIOLOGY

PNAS Plus: Defects in coatamer protein I (COPI) transport cause blood feeding-induced mortality in Yellow Fever mosquitoes

Jun Isoe, Jennifer Collins, Hemant Badgandi, W. Anthony Day, and Roger L. Miesfeld
PNAS, Jun 2011; 108: E211 - E217.

©2011 LI-COR, Inc. LI-COR, Odyssey, MousePOD, IRDye, and In-Cell Western are trademarks and registered trademarks of LI-COR, Inc., in the United States and other countries. All other trademarks belong to their respective owners.



4647 Superior St. • P.O. Box 4000 • Lincoln, Nebraska 68504
North America: 800-645-4267 • International: 402-467-0700

LI-COR GmbH, Germany, Serving Europe, Middle East and Africa: +49 (0) 6172 17 17 771

LI-COR Ltd, UK, Serving UK, Ireland and Scandinavia: +44 (0) 1223 422104

All other countries, contact LI-COR Biosciences or a local LI-COR distributor:
<http://www.licor.com/distributors>

www.licor.com/bio

Doc #979-07647
Rev H 0811