

Odyssey® Publications

Selected from 242 publications in July 1, 2011 – September 30, 2011, HighWire

Fall 2011

Volume 13

TABLE OF CONTENTS

	Page
Apoptosis	2
Autophagy	2
Angiogenesis	2
Cancer	2
Cardiovascular	4
Cell Adhesion, Migration and Invasion	4
Cell Cycle/DNA Replication	5
Cell Signaling	5
Diabetes/Metabolism/Obesity	6
Immunology	7
Ion Channels and Transporters	8
Micro RNA	9
Muscle	9
Neuroscience	10
Nephrology/Kidney	10
Phosphatases	10
Proteases	11
Receptors	11
RNAi	12
ROS – Reactive Oxygen Species	13
Transcription Factors	13
Ubiquitination and Sumoylation	14
Virology/Microbiology	14
Applications	16
<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® Accessory</i>	
<i>Reverse Phase Protein Arrays</i>	
<i>RNA/Protein Interactions</i>	
<i>siRNA Screening</i>	
<i>Western (Quantitative, Molecules/Cell)</i>	
PNAS Publications This Quarter	18

LI-COR®

Experience Excellence

PUBLICATIONS REFERENCING ODYSSEY DATA

RESEARCH AREAS

APOPTOSIS

Targeted Deletion of Adipocytes by Apoptosis Leads to Adipose Tissue Recruitment of Alternatively Activated M2 Macrophages

Pamela Fischer-Posovszky, Qiong A. Wang, Ingrid Wernstedt Asterholm, Joseph M. Rutkowski, and Philipp E. Scherer
Endocrinology, Aug 2011; 152: 3074 - 3081.

Novel Mechanism of Anti-apoptotic Function of 78-kDa Glucose-regulated Protein (GRP78): ENDOCRINE RESISTANCE FACTOR IN BREAST CANCER, THROUGH RELEASE OF B-CELL LYMPHOMA 2 (BCL-2) FROM BCL-2-INTERACTING KILLER (BIK)

Hui Zhou, Yi Zhang, Yong Fu, Lauren Chan, and Amy S. Lee
J. Biol. Chem., Jul 2011; 286: 25687 - 25696.

Role for hACF1 in the G2/M damage checkpoint

Sara Sánchez-Molina, Oliver Mortusewicz, Béatrice Bieber, Susanne Auer, Maren Eckey, Heinrich Leonhardt, Anna A. Friedl, and Peter B. Becker

Nucleic Acids Res., Jul 2011; 10.1093/nar/gkr435.

IRES-mediated translation of cellular messenger RNA operates in eIF2 α - independent manner during stress

Nehal Thakor and Martin Holcik

Nucleic Acids Res., Sep 2011; 10.1093/nar/gkr701.

AUTOPHAGY

An Atg13 Protein-mediated Self-association of the Atg1 Protein Kinase Is Important for the Induction of Autophagy

Yuh-Ying Yeh, Khyati H. Shah, and Paul K. Herman

J. Biol. Chem., Aug 2011; 286: 28931 - 28939.

ANGIOGENESIS

Endorepellin, the Angiostatic Module of Perlecan, Interacts with Both the $\alpha 2\beta 1$ Integrin and Vascular Endothelial Growth Factor Receptor 2 (VEGFR2): A DUAL RECEPTOR ANTAGONISM

Atul Goyal, Nutan Pal, Matthew Concannon, Matthew Paul, Mike Doran, Chiara Poluzzi, Kiyotoshi Sekiguchi, John M. Whitelock, Thomas Neill, and Renato V. Iozzo

J. Biol. Chem., Jul 2011; 286: 25947 - 25962.

CANCER

Epigenetic Regulation of Phosphatidylinositol 3,4,5-Triphosphate-dependent Rac Exchanger 1 Gene Expression in Prostate Cancer Cells

Chuu-Yun A. Wong, Hada Wuriyanghan, Yan Xie, Ming-Fong Lin, Peter W. Abel, and Yaping Tu

J. Biol. Chem., Jul 2011; 286: 25813 - 25822.

Mechanisms of cytotoxicity to Pim kinase inhibitor, SGI-1776, in acute myeloid leukemia

Lisa S. Chen, Sanjeev Redkar, Pietro Taverna, Jorge E. Cortes, and Varsha Gandhi

Blood, Jul 2011; 118: 693 - 702.

Rapamycin Partially Mimics the Anticancer Effects of Calorie Restriction in a Murine Model of Pancreatic Cancer

Laura M. Lashinger, Lauren M. Malone, Graham W. Brown, Elizabeth A. Daniels, Jason A. Goldberg, Glen Otto, Susan M. Fischer, and Stephen D. Hursting

Cancer Prevention Research, Jul 2011; 4: 1041 - 1051.

Genetic Reduction of Insulin-like Growth Factor-1 Mimics the Anticancer Effects of Calorie Restriction on Cyclooxygenase-2–Driven Pancreatic Neoplasia

Laura M. Lashinger, Lauren M. Malone, Mark J. McArthur, Jason A. Goldberg, Elizabeth A. Daniels, Amy Pavone, Jennifer K. Colby, Nicole C. Smith, Susan N. Perkins, Susan M. Fischer, and Stephen D. Hursting
Cancer Prevention Research, Jul 2011; 4: 1030 - 1040.

Killing of Chronic Lymphocytic Leukemia by the Combination of Fludarabine and Oxaliplatin Is Dependent on the Activity of XPF Endonuclease

Alma Zecevic, Deepa Sampath, Brett Ewald, Rong Chen, William Wierda, and William Plunkett
Clin. Cancer Res., Jul 2011; 17: 4731 - 4741.

Truncated ErbB2 Expressed in Tumor Cell Nuclei Contributes to Acquired Therapeutic Resistance to ErbB2 Kinase Inhibitors

Wenle Xia, Zuguo Liu, Rongrong Zong, Leihua Liu, Sumin Zhao, Sarah S. Bacus, Yubin Mao, Jia He, Julia D. Wulfkuhle, Emanuel F. Petricoin, III, Takuya Osada, Xiao-Yi Yang, Zachary C. Hartman, Timothy M. Clay, Kimberly L. Blackwell, Herbert K. Lyerly, and Neil L. Spector
Mol. Cancer Ther., Aug 2011; 10: 1367 - 1374.

Thioredoxin-Interacting Protein Mediates TRX1 Translocation to the Plasma Membrane in Response to Tumor Necrosis Factor- α : A Key Mechanism for Vascular Endothelial Growth Factor Receptor-2 Transactivation by Reactive Oxygen Species

Cameron World, Oded N. Spindel, and Bradford C. Berk
Arterioscler Thromb Vasc Biol, Aug 2011; 31: 1890 - 1897.

Peloruside- and Laulimalide-Resistant Human Ovarian Carcinoma Cells Have β I-Tubulin Mutations and Altered Expression of β II- and β III-Tubulin Isoforms

Arun Kanakkanthara, Anja Wilmes, Aurora O'Brate, Daniel Escuin, Ariane Chan, Ada Gjyzezi, Janet Crawford, Pisana Rawson, Bronwyn Kivell, Peter T. Northcote, Ernest Hamel, Paraskevi Giannakakou, and John H. Miller
Mol. Cancer Ther., Aug 2011; 10: 1419 - 1429.

Berberine Suppresses Androgen Receptor Signaling in Prostate Cancer

Jing Li, Bo Cao, Xichun Liu, Xueqi Fu, Zhenggang Xiong, Li Chen, Oliver Sartor, Yan Dong, and Haitao Zhang
Mol. Cancer Ther., Aug 2011; 10: 1346 - 1356.

A *Cyp2a* polymorphism predicts susceptibility to NNK-induced lung tumorigenesis in mice

M.Christine Hollander, Xin Zhou, Colleen R. Maier, Andrew D. Patterson, Xinxin Ding, and Phillip A. Dennis
Carcinogenesis, Aug 2011; 32: 1279 - 1284.

Engineered Bivalent Ligands to Bias ErbB Receptor-mediated Signaling and Phenotypes

Steven M. Jay, Elma Kurtagic, Luis M. Alvarez, Seymour de Picciotto, Edgar Sanchez, Jessica F. Hawkins, Robin N. Prince, Yadir Guerrero, Carolyn L. Treasure, Richard T. Lee, and Linda G. Griffith
J. Biol. Chem., Aug 2011; 286: 27729 - 27740.

Inhibition of gap junction transfer sensitizes thyroid cancer cells to anoikis

Kirk Jensen, Aneeta Patel, Joanna Klubo-Gwiezdzinska, Andrew Bauer, and Vasyl Vasko
Endocr. Relat. Cancer, Sep 2011; 18: 613 - 626.

Activating Enhancer-Binding Protein-2 β Nucleolar Localization Predicts Poor Survival After Stage I Non-Small Cell Lung Cancer Resection

Min P. Kim, Ying Chen, B. Nebiyu Bekele, Adriana Lopez, Abha Khanna, Jie Qing Chen, Margaret R. Spitz, Carmen Behrens, Luisa Solis, Marnie Wismach, Lin Ji, Ignacio I. Wistuba, Jack A. Roth, and Ruth L. Katz
Ann. Thorac. Surg., Sep 2011; 92: 1044 – 1050.

Insulin Increases *De Novo* Steroidogenesis in Prostate Cancer Cells

Amy A. Lubik, Jennifer H. Gunter, Stephen C. Hendy, Jennifer A. Locke, Hans H. Adomat, Vanessa Thompson, Adrian Herington, Martin E. Gleave, Michael Pollak, and Colleen C. Nelson
Cancer Res., Sep 2011; 71: 5754 - 5764.

Inhibition of gap junction transfer sensitizes thyroid cancer cells to anoikis

Kirk Jensen, Aneeta Patel, Joanna Klubo-Gwiezdzinska, Andrew Bauer, and Vasyl Vasko
Endocr. Relat. Cancer, Sep 2011; 18: 613 - 626.

Potent, Selective Inhibitors of Fibroblast Growth Factor Receptor Define Fibroblast Growth Factor Dependence in Preclinical Cancer Models

Matthew Squires, George Ward, Gordan Saxty, Valerio Berdini, Anne Cleasby, Peter King, Patrick Angibaud, Tim Perera, Lynsey Fazal, Douglas Ross, Charlotte Griffiths Jones, Andrew Madin, Rajdeep K. Benning, Emma Vickerstaffe, Alistair O'Brien, Martyn Frederickson, Michael Reader, Christopher Hamlett, Michael A. Batey, Sharna Rich, Maria Carr, Darcey Miller, Ruth Feltell, Abarna Thiru, Susanne Bethell, Lindsay A. Devine, Brent L. Graham, Andrew Pike, Jose Cosme, Edward J. Lewis, Eddy Freyne, John Lyons, Julie Irving, Christopher Murray, David R. Newell, and Neil T. Thompson
Mol. Cancer Ther., Sep 2011; 10: 1542 - 1552.

Clusterin Inhibition Using OGX-011 Synergistically Enhances Hsp90 Inhibitor Activity by Suppressing the Heat Shock Response in Castrate-Resistant Prostate Cancer

Francois Lamoureux, Christian Thomas, Min-Jean Yin, Hidetoshi Kuruma, Eliana Beraldi, Ladan Fazli, Amina Zoubeidi, and Martin E. Gleave
Cancer Res., Sep 2011; 71: 5838 - 5849.

Novel Positron Emission Tomography Tracer Distinguishes Normal from Cancerous Cells

Muhammad Saeed, David Sheff, and Amnon Kohen
J. Biol. Chem., Sep 2011; 286: 33872 - 33878.

Taxane-Induced Blockade to Nuclear Accumulation of the Androgen Receptor Predicts Clinical Responses in Metastatic Prostate Cancer

Medha S. Darshan, Matthew S. Loftus, Maria Thadani-Mulero, Benjamin P. Levy, Daniel Escuin, Xi Kathy Zhou, Ada Gjyzezi, Chantal Chanel-Vos, Ruoqian Shen, Scott T. Tagawa, Neil H. Bander, David M. Nanus, and Paraskevi Giannakakou
Cancer Res., Sep 2011; 71: 6019 - 6029.

Cardiovascular

Doxycycline, a matrix metalloprotease inhibitor, reduces vascular remodeling and damage after cerebral ischemia in stroke-prone spontaneously hypertensive rats

Paulo W. Pires, Curt T. Rogers, Jonathon L. McClain, Hannah S. Garver, Gregory D. Fink, and Anne M. Dorrance
Am J Physiol Heart Circ Physiol, Jul 2011; 301: H87 - H97.

Monocytic microparticles activate endothelial cells in an IL-1 β -dependent manner

Jian-Guo Wang, Julie C. Williams, Beckley K. Davis, Ken Jacobson, Claire M. Doerschuk, Jenny P.-Y. Ting, and Nigel Mackman
Blood, Aug 2011; 118: 2366 - 2374.

Contributions of VEGF to age-dependent transmural gradients in contractile protein expression in ovine carotid arteries

Stacy M. Butler, Jenna M. Abrassart, Margaret C. Hubbell, Olayemi Adeoye, Andrew Semotiuk, James M. Williams, Eugenia Mata-Greenwood, Omid Khorram, and William J. Pearce
Am J Physiol Cell Physiol, Sep 2011; 301: C653 - C666.

Cell Adhesion, Migration and Invasion

Exocyst Complex Component 3-like 2 (EXOC3L2) Associates with the Exocyst Complex and Mediates Directional Migration of Endothelial Cells

Irmeli Barkefors, Peder Fredlund Fuchs, Johan Heldin, Tobias Bergström, Karin Forsberg-Nilsson, and Johan Kreuger
J. Biol. Chem., Jul 2011; 286: 24189 - 24199.

CD82 gene suppression in endometrial stromal cells leads to increase of the cell invasiveness in the endometriotic milieu

Ming-Qing Li, Xiao-Fan Hou, Shi-Jian Lv, Yu-Han Meng, Xiao-Qiu Wang, Chuan-Ling Tang, and Da-Jin Li
J. Mol. Endocrinol., Aug 2011; 47: 195 - 208.

Hepatitis C Virus Infection Causes Cell Cycle Arrest at the Level of Initiation of Mitosis

Rathi P. Kannan, Lucinda L. Hensley, Lauren E. Evers, Stanley M. Lemon, and David R. McGivern
J. Virol., Aug 2011; 85: 7989 - 8001.

Proteolytic Activation of Pro-Macrophage-Stimulating Protein by Hepsin

Rajkumar Ganesan, Ganesh A. Kolumam, S. Jack Lin, Ming-Hong Xie, Lydia Santell, Thomas D. Wu, Robert A. Lazarus, Amitabha Chaudhuri, and Daniel Kirchhofer
Mol. Cancer Res., Sep 2011; 9: 1175 - 1186.

Cell Cycle/DNA Replication**Role for hACF1 in the G2/M damage checkpoint**

Sara Sánchez-Molina, Oliver Mortusewicz, Béatrice Bieber, Susanne Auer, Maren Eckey, Heinrich Leonhardt, Anna A. Friedl, and Peter B. Becker
Nucleic Acids Res., Jul 2011; 10.1093/nar/gkr435.

Hepatitis C Virus Infection Causes Cell Cycle Arrest at the Level of Initiation of Mitosis

Rathi P. Kannan, Lucinda L. Hensley, Lauren E. Evers, Stanley M. Lemon, and David R. McGivern
J. Virol., Aug 2011; 85: 7989 - 8001.

Overlapping kinetochore targets of CK2 and Aurora B kinases in mitotic regulation

Yutian Peng, Catherine C. L. Wong, Yuko Nakajima, Randall G. Tyers, Ali S. Sarkeshik, John Yates, III, David G. Drubin, and Georjana Barnes
Mol. Biol. Cell, Aug 2011; 22: 2680 - 2689.

Time-Dependent Quantitative Multicomponent Control of the G₁-S Network by the Stress-Activated Protein Kinase Hog1 upon Osmostress

Miquel Àngel Adrover, Zhike Zi, Alba Duch, Jörg Schaber, Alberto González-Novo, Javier Jimenez, Mariona Nadal-Ribelles, Josep Clotet, Edda Klipp, and Francesc Posas
Sci. Signal., Sep 2011; 4: ra63.

Histone hypoacetylation is required to maintain late replication timing of constitutive heterochromatin

Corella S. Casas-Delucchi, Joke G. van Bommel, Sebastian Haase, Henry D. Herce, Danny Nowak, Daniela Meilinger, Jeffrey H. Stear, Heinrich Leonhardt, and M. Cristina Cardoso
Nucleic Acids Res., Sep 2011; 10.1093/nar/gkr723.

Cell Signaling**TREM-1 Amplifies Corneal Inflammation after *Pseudomonas aeruginosa* Infection by Modulating Toll-Like Receptor Signaling and Th1/Th2-Type Immune Responses**

Minhao Wu, Anping Peng, Mingxia Sun, Qiuchan Deng, Linda D. Hazlett, Jin Yuan, Xialin Liu, Qianying Gao, Lianqiang Feng, Junfang He, Ping Zhang, and Xi Huang
Infect. Immun., Jul 2011; 79: 2709 - 2716.

Involvement of Epidermal Growth Factor Receptor Signaling in Estrogen Inhibition of Oocyte Maturation Mediated Through the G Protein-Coupled Estrogen Receptor (Gper) in Zebrafish (*Danio rerio*)

Candace Peyton and Peter Thomas
Biol. Reprod., Jul 2011; 85: 42 - 50.

From Molecular Details of the Interplay between Transmembrane Helices of the Thyrotropin Receptor to General Aspects of Signal Transduction in Family A G-protein-coupled Receptors (GPCRs)

Gunnar Kleinau, Inna Hoyer, Annika Kreuchwig, Ann-Karin Haas, Claudia Rutz, Jens Ferkert, Catherine L. Worth, Gerd Krause, and Ralf Schüle
J. Biol. Chem., Jul 2011; 286: 25859 – 25871.

The PI3K/Akt pathway is involved in early infection of some exogenous avian leukosis viruses

Shao-zhen Feng, Wei-sheng Cao, and Ming Liao
J. Gen. Virol., Jul 2011; 92: 1688 - 1697.

Ca²⁺/Calmodulin-dependent Protein Kinase Kinase β Is Regulated by Multisite Phosphorylation

Michelle F. Green, John W. Scott, Rohan Steel, Jonathan S. Oakhill, Bruce E. Kemp, and Anthony R. Means
J. Biol. Chem., Aug 2011; 286: 28066 - 28079.

Processive phosphorylation of ERK MAP kinase in mammalian cells

Kazuhiro Aoki, Masashi Yamada, Katsuyuki Kunida, Shuhei Yasuda, and Michiyuki Matsuda
PNAS, Aug 2011; 108: 12675 - 12680.

Engineered Bivalent Ligands to Bias ErbB Receptor-mediated Signaling and Phenotypes

Steven M. Jay, Elma Kurtagic, Luis M. Alvarez, Seymour de Picciotto, Edgar Sanchez, Jessica F. Hawkins, Robin N. Prince, Yadir Guerrero, Carolyn L. Treasure, Richard T. Lee, and Linda G. Griffith
J. Biol. Chem., Aug 2011; 286: 27729 - 27740.

TGF- β /Smad3 Signaling Promotes Renal Fibrosis by Inhibiting miR-29

Wei Qin, Arthur C.K. Chung, Xiao R. Huang, Xiao-Ming Meng, David S.C. Hui, Cheuk-Man Yu, Joseph J.Y. Sung, and Hui Y. Lan
J. Am. Soc. Nephrol., Aug 2011; 22: 1462 - 1474.

An Atg13 Protein-mediated Self-association of the Atg1 Protein Kinase Is Important for the Induction of Autophagy

Yuh-Ying Yeh, Khyati H. Shah, and Paul K. Herman
J. Biol. Chem., Aug 2011; 286: 28931 - 28939.

NMDA Receptor Signaling in Oligodendrocyte Progenitors Is Not Required for Oligodendrogenesis and Myelination

Lindsay M. De Biase, Shin H. Kang, Emily G. Baxi, Masahiro Fukaya, Michele L. Pucak, Masayoshi Mishina, Peter A. Calabresi, and Dwight E. Bergles
J. Neurosci., Aug 2011; 31: 12650 - 12662.

Src Homology Domain 2-containing Protein-tyrosine Phosphatase-1 (SHP-1) Binds and Dephosphorylates G α -interacting, Vesicle-associated Protein (GIV)/Girdin and Attenuates the GIV-Phosphatidylinositol 3-Kinase (PI3K)-Akt Signaling Pathway

Yash Mittal, Yelena Pavlova, Mikel Garcia-Marcos, and Pradipta Ghosh
J. Biol. Chem., Sep 2011; 286: 32404 - 32415.

β -catenin tyrosine 654 phosphorylation increases Wnt signalling and intestinal tumorigenesis

Wendy van Veelen, Ngoc H Le, Werner Helvensteijn, Lau Blonden, Myrte Theeuwes, Elvira R M Bakker, Patrick F Franken, Léon van Gurp, Frits Meijlink, Martin A van der Valk, Ernst J Kuipers, Riccardo Fodde, and Ron Smits
Gut, Sep 2011; 60: 1204 - 1212.

Gender-Specific Interplay of Signaling through β -Catenin and CAR in the Regulation of Xenobiotic-Induced Hepatocyte Proliferation

Albert Braeuning, Yvonne Heubach, Thomas Knorpp, Marta Anna Kowalik, Markus Templin, Amedeo Columbano, and Michael Schwarz
Toxicol. Sci., Sep 2011; 123: 113 - 122.

WNT Protein-independent Constitutive Nuclear Localization of β -Catenin Protein and Its Low Degradation Rate in Thalamic Neurons

Katarzyna Misztal, Marta B. Wisniewska, Mateusz Ambrozkiwicz, Andrzej Nagalski, and Jacek Kuznicki
J. Biol. Chem., Sep 2011; 286: 31781 - 31788.

Diabetes/ Metabolism/Obesity

Ovine corpus luteum proteins, with functions including oxidative stress and lipid metabolism, show complex alterations during implantation

Mitra Arianmanesh, Rebecca H McIntosh, Richard G Lea, Paul A Fowler, and Kaïs H Al-Gubory
J. Endocrinol., Jul 2011; 210: 47 - 58.

Insulinotropic and Antidiabetic Effects of 17 β -Estradiol and the GPR30 Agonist G-1 on Human Pancreatic Islets

Rajesh Kumar, Alexander Balhuizen, Stefan Amisten, Ingmar Lundquist, and Albert Salehi
Endocrinology, Jul 2011; 152: 2568 - 2579.

Klotho Depletion Contributes to Increased Inflammation in Kidney of the *db/db* Mouse Model of Diabetes via RelA (Serine)⁵³⁶ Phosphorylation

Yanhua Zhao, Srijita Banerjee, Nilay Dey, Wanda S. LeJeune, Partha S. Sarkar, Reynolds Brobey, Kevin P. Rosenblatt, Ronald G. Tilton, and Sanjeev Choudhary
Diabetes, Jul 2011; 60: 1907 - 1916.

Chemerin, a Novel Peroxisome Proliferator-activated Receptor γ (PPAR γ) Target Gene That Promotes Mesenchymal Stem Cell Adipogenesis

Shanmugam Muruganandan, Sebastian D. Parlee, Jillian L. Rourke, Matthew C. Ernst, Kerry B. Goralski, and Christopher J. Sinal
J. Biol. Chem., Jul 2011; 286: 23982 - 23995.

Maternal Obesity-Impaired Insulin Signaling in Sheep and Induced Lipid Accumulation and Fibrosis in Skeletal Muscle of Offspring

Xu Yan, Yan Huang, Jun-Xing Zhao, Nathan M. Long, Adam B. Uthlaut, Mei-Jun Zhu, Stephen P. Ford, Peter W. Nathanielsz, and Min Du
Biol. Reprod., Jul 2011; 85: 172 - 178.

A lipid-droplet-targeted O-GlcNAcase isoform is a key regulator of the proteasome

Chithra N. Keembiyehetty, Anna Krzeslak, Dona C. Love, and John A. Hanover
J. Cell Sci., Aug 2011; 124: 2851 - 2860.

Folate Intake, *Mthfr* Genotype, and Sex Modulate Choline Metabolism in Mice

Tina W. Chew, Xinyin Jiang, Jian Yan, Wei Wang, Amanda L. Lusa, Bradley J. Carrier, Allyson A. West, Olga V. Malysheva, J. Thomas Brenna, Jesse F. Gregory, III, and Marie A. Caudill
J. Nutr., Aug 2011; 141: 1475 - 1481.

Targeted Deletion of Adipocytes by Apoptosis Leads to Adipose Tissue Recruitment of Alternatively Activated M2 Macrophages

Pamela Fischer-Posovszky, Qiong A. Wang, Ingrid Wernstedt Asterholm, Joseph M. Rutkowski, and Philipp E. Scherer
Endocrinology, Aug 2011; 152: 3074 - 3081.

Akt-Dependent Glucose Metabolism Promotes Mcl-1 Synthesis to Maintain Cell Survival and Resistance to Bcl-2 Inhibition

Jonathan L. Coloff, Andrew N. Macintyre, Amanda G. Nichols, Tingyu Liu, Catherine A. Gallo, David R. Plas, and Jeffrey C. Rathmell
Cancer Res., Aug 2011; 71: 5204 - 5213.

Functional Plasticity of Central TRPV1 Receptors in Brainstem Dorsal Vagal Complex Circuits of Streptozotocin-Treated Hyperglycemic Mice

Andrea Zsombok, Muthu D. Bhaskaran, Hong Gao, Andrei V. Derbenev, and Bret N. Smith
J. Neurosci., Sep 2011; 31: 14024 - 14031.

The Unconventional Role of Acid Sphingomyelinase in Regulation of Retinal Microangiopathy in Diabetic Human and Animal Models

Madalina Opreanu, Maria Tikhonenko, Svetlana Bozack, Todd A. Lydic, Gavin E. Reid, Kelly M. McSorley, Andrew Sochacki, Gloria I. Perez, Walter J. Esselman, Timothy Kern, Richard Kolesnick, Maria B. Grant, and Julia V. Busik
Diabetes, Sep 2011; 60: 2370 - 2378.

Immunology

TREM-1 Amplifies Corneal Inflammation after *Pseudomonas aeruginosa* Infection by Modulating Toll-Like Receptor Signaling and Th1/Th2-Type Immune Responses

Minhao Wu, Anping Peng, Mingxia Sun, Qiuchan Deng, Linda D. Hazlett, Jin Yuan, Xialin Liu, Qianying Gao, Lianqiang Feng, Junfang He, Ping Zhang, and Xi Huang
Infect. Immun., Jul 2011; 79: 2709 - 2716.

Retinoic Acid-induced Gene-I (RIG-I) Associates with Nucleotide-binding Oligomerization Domain-2 (NOD2) to Negatively Regulate Inflammatory Signaling

Stefanie A. Morosky, Jianzhong Zhu, Amitava Mukherjee, Saumendra N. Sarkar, and Carolyn B. Coyne
J. Biol. Chem., Aug 2011; 286: 28574 - 28583.

Molecular Mechanism of Signal Perception and Integration by the Innate Immune Sensor Retinoic Acid-inducible Gene-I (RIG-I)

Marco Binder, Florian Eberle, Stefan Seitz, Norbert Mücke, Christian M. Hüber, Narsis Kiani, Lars Kaderali, Volker Lohmann, Alexander Dalpke, and Ralf Bartenschlager
J. Biol. Chem., Aug 2011; 286: 27278 - 27287.

Ion Channels and Transporters

A haploid genetic screen identifies the major facilitator domain containing 2A (MFSD2A) transporter as a key mediator in the response to tunicamycin

Jan H. Reiling, Clary B. Clish, Jan E. Carette, Malini Varadarajan, Thijn R. Brummelkamp, and David M. Sabatini
PNAS, Jul 2011; 108: 11756 - 11765.

Probing Conformational Rescue Induced by a Chemical Corrector of F508del-Cystic Fibrosis Transmembrane Conductance Regulator (CFTR) Mutant

Wilson Yu, Patrick Kim Chiaw, and Christine E. Bear
J. Biol. Chem., Jul 2011; 286: 24714 - 24725.

Placental Expression of the Heme Transporter, Feline Leukemia Virus Subgroup C Receptor, Is related to Maternal Iron Status in Pregnant Adolescents

Lindsay M. Jaacks, Melissa F. Young, Bridget V. Essley, Thomas J. McNanley, Elizabeth M. Cooper, Eva K. Pressman, Allison W. McIntyre, Mark S. Orlando, Janis L. Abkowitz, Ronnie Guillet, and Kimberly O. O'Brien
J. Nutr., Jul 2011; 141: 1267 - 1272.

Amino acid absorption and homeostasis in mice lacking the intestinal peptide transporter PEPT1

Anna-Maria Nässl, Isabel Rubio-Aliaga, Henning Fenselau, Mena Katharina Marth, Gabor Kottra, and Hannelore Daniel
Am J Physiol Gastrointest Liver Physiol, Jul 2011; 301: G128 - G137.

Trafficking-deficient hERG K⁺ channels linked to long QT syndrome are regulated by a microtubule-dependent quality control compartment in the ER

Jennifer L. Smith, Christie M. McBride, Parvathi S. Nataraj, Daniel C. Bartos, Craig T. January, and Brian P. Delisle
Am J Physiol Cell Physiol, Jul 2011; 301: C75 - C85.

Chloride Channel (Clc)-5 Is Necessary for Exocytic Trafficking of Na⁺/H⁺ Exchanger 3 (NHE3)

Zhihong Lin, Shi Jin, Xiaohong Duan, Tong Wang, Sabrina Martini, Phuson Hulamm, Boyoung Cha, Ann Hubbard, Mark Donowitz, and Sandra E. Guggino
J. Biol. Chem., Jul 2011; 286: 22833 - 22845.

Transmembrane Helix 1 Contributes to Substrate Translocation and Protein Stability of Bile Acid Transporter SLC10A2

Tatiana Claro da Silva, Naissan Hussainzada, Chandra M. Khantwal, James E. Polli, and Peter W. Swaan
J. Biol. Chem., Aug 2011; 286: 27322 - 27332.

VEGF regulates TRPC6 channels in podocytes

Florian Thilo, Ying Liu, Christoph Loddenkemper, Ralf Schuelein, Antje Schmidt, Zengcheng Yan, Zhiming Zhu, Andreas Zakrzewicz, Maik Gollasch, and Martin Tepel
Nephrol. Dial. Transplant., Aug 2011; 10.1093/ndt/gfr457.

Erythropoietin Increases Expression and Function of Transient Receptor Potential Canonical 5 Channels

Ying Liu, Yunfei Xu, Florian Thilo, Ulla G. Friis, Boye L. Jensen, Alexandra Scholze, Junhua Zheng, and Martin Tepel
Hypertension, Aug 2011; 58: 317 - 324.

Stanniocalcin 2 Is a Negative Modulator of Store-Operated Calcium Entry

William Zeiger, Daisuke Ito, Carol Swetlik, Masatsugu Oh-hora, Mitchel L. Villereal, and Gopal Thinakaran
Mol. Cell. Biol., Sep 2011; 31: 3710 - 3722.

Micro RNA

TGF- β /Smad3 Signaling Promotes Renal Fibrosis by Inhibiting miR-29

Wei Qin, Arthur C.K. Chung, Xiao R. Huang, Xiao-Ming Meng, David S.C. Hui, Cheuk-Man Yu, Joseph J.Y. Sung, and Hui Y. Lan
J. Am. Soc. Nephrol., Aug 2011; 22: 1462 - 1474.

The DExD/H box ATPase Dhh1 functions in translational repression, mRNA decay, and processing body dynamics

Johanna S. Carroll, Sarah E. Munchel, and Karsten Weis
J. Cell Biol., Aug 2011; 194: 527 - 537.

Muscle

Muscle pathology without severe nerve pathology in a new mouse model of Charcot–Marie–Tooth disease type 2E

Hailian Shen, Devin M. Barry, Jeffrey M. Dale, Virginia B. Garcia, Nigel A. Calcutt, and Michael L. Garcia
Hum. Mol. Genet., Jul 2011; 20: 2535 - 2548.

Maternal Obesity-Impaired Insulin Signaling in Sheep and Induced Lipid Accumulation and Fibrosis in Skeletal Muscle of Offspring

Xu Yan, Yan Huang, Jun-Xing Zhao, Nathan M. Long, Adam B. Uthlaut, Mei-Jun Zhu, Stephen P. Ford, Peter W. Nathanielsz, and Min Du
Biol Reprod, Jul 2011; 85: 172 - 178.

Transgenic overexpression of the $\alpha 7$ integrin reduces muscle pathology and improves viability in the dy^w mouse model of merosin-deficient congenital muscular dystrophy type 1A

Jinger A. Doe, Ryan D. Wuebbles, Erika T. Allred, Jachinta E. Rooney, Margaret Elorza, and Dean J. Burkin
J. Cell Sci., Jul 2011; 124: 2287 - 2297.

Reversible molecular pathology of skeletal muscle in spinal muscular atrophy

Chantal A. Mutsaers, Thomas M. Wishart, Douglas J. Lamont, Markus Riessland, Julia Schreml, Laura H. Comley, Lyndsay M. Murray, Simon H. Parson, Hanns Lochmüller, Brunhilde Wirth, Kevin Talbot, and Thomas H. Gillingwater
Hum. Mol. Genet., Aug 2011; 10.1093/hmg/ddr360.

COOH-terminal truncation of flightin decreases myofilament lattice organization, cross-bridge binding, and power output in *Drosophila* indirect flight muscle

Bertrand C. W. Tanner, Mark S. Miller, Becky M. Miller, Panagiotis Lekkas, Thomas C. Irving, David W. Maughan, and Jim O. Vigoreaux
Am J Physiol Cell Physiol, Aug 2011; 301: C383 - C391.

Immunocytochemical evidence for PDBu-induced activation of RhoA/ROCK in human internal anal sphincter smooth muscle cells

Jagmohan Singh, Pinckney J. Maxwell, IV, and Satish Rattan
Am J Physiol Gastrointest Liver Physiol, Aug 2011; 301: G317 - G325.

Vegetarians have a reduced skeletal muscle carnitine transport capacity

Francis B Stephens, Kanagaraj Marimuthu, Yi Cheng, Nitin Patel, Despina Constantin, Elizabeth J Simpson, and Paul L Greenhaff
Am J Clin Nutr, Sep 2011; 94: 938 – 944.

Effective fiber hypertrophy in satellite cell-depleted skeletal muscle

John J. McCarthy, Jyothi Mula, Mitsunori Miyazaki, Rod Erfani, Kelcye Garrison, Amreen B. Farooqui, Ratchakrit Srikuea, Benjamin A. Lawson, Barry Grimes, Charles Keller, Gary Van Zant, Kenneth S. Campbell, Karyn A. Esser, Esther E. Dupont-Versteegden, and Charlotte A. Peterson
Development, Sep 2011; 138: 3657 - 3666.

Neuroscience

CNS expression of glucocerebrosidase corrects α -synuclein pathology and memory in a mouse model of Gaucher-related synucleinopathy

S. Pablo Sardi, Jennifer Clarke, Cathrine Kinnecom, Thomas J. Tamsett, Lingyun Li, Lisa M. Stanek, Marco A. Passini, Gregory A. Grabowski, Michael G. Schlossmacher, Richard L. Sidman, Seng H. Cheng, and Lamya S. Shihabuddin
PNAS, Jul 2011; 108: 12101 - 12106.

S-Nitrosylation activates Cdk5 and contributes to synaptic spine loss induced by β -amyloid peptide

Jing Qu, Tomohiro Nakamura, Gang Cao, Emily A. Holland, Scott R. McKercher, and Stuart A. Lipton
PNAS, Aug 2011; 108: 14330 - 14335.

Blast-induced phenotypic switching in cerebral vasospasm

Patrick W. Alford, Borna E. Dabiri, Josue A. Goss, Matthew A. Hemphill, Mark D. Brigham, and Kevin Kit Parker
PNAS, Aug 2011; 108: 12705 - 12710.

Permanent deficits in brain functions caused by long-term ketamine treatment in mice

L Sun, WP Lam, YW Wong, LH Lam, HC Tang, MS Wai, YT Mak, F Pan, and DTYew
Human and Experimental Toxicology, Sep 2011; 30: 1287 - 1296.

Nephrology/Kidney

Distinct Actions of Endothelin A-Selective Versus Combined Endothelin A/B Receptor Antagonists in Early Diabetic Kidney Disease

Mohamed A. Saleh, Jennifer S. Pollock, and David M. Pollock
J. Pharmacol. Exp. Ther., Jul 2011; 338: 263 - 270.

VEGF regulates TRPC6 channels in podocytes

Florian Thilo, Ying Liu, Christoph Loddenkemper, Ralf Schuelein, Antje Schmidt, Zengcheng Yan, Zhiming Zhu, Andreas Zakrzewicz, Maik Gollasch, and Martin Tepel
Nephrol. Dial. Transplant., Aug 2011; 10.1093/ndt/gfr457.

TGF- β /Smad3 Signaling Promotes Renal Fibrosis by Inhibiting miR-29

Wei Qin, Arthur C.K. Chung, Xiao R. Huang, Xiao-Ming Meng, David S.C. Hui, Cheuk-Man Yu, Joseph J.Y. Sung, and Hui Y. Lan
J. Am. Soc. Nephrol., Aug 2011; 22: 1462 - 1474.

Renal inflammation and elevated blood pressure in a mouse model of reduced β -ENaC

Heather A. Drummond, Samira C. Grifoni, Ahmed Abu-Zaid, Monette Gousset, Rumbidayi Chiposi, John M. Barnard, Beau Murphey, and David E. Stec
Am J Physiol Renal Physiol, Aug 2011; 301: F443 - F449.

Functional Plasticity of Central TRPV1 Receptors in Brainstem Dorsal Vagal Complex Circuits of Streptozotocin-Treated Hyperglycemic Mice

Andrea Zsombok, Muthu D. Bhaskaran, Hong Gao, Andrei V. Derbenev, and Bret N. Smith
J. Neurosci., Sep 2011; 31: 14024 - 14031.

Phosphatases

Microarray and Proteomic Analysis of Breast Cancer Cell and Osteoblast Co-cultures: ROLE OF OSTEOBLAST MATRIX METALLOPROTEINASE (MMP)-13 IN BONE METASTASIS

Charlotte Morrison, Stephanie Mancini, Jane Cipollone, Reinhild Kappelhoff, Calvin Roskelley, and Christopher Overall
J. Biol. Chem., Sep 2011; 286: 34271 - 34285.

Inhibition of Protein-tyrosine Phosphatase 1B (PTP1B) Mediates Ubiquitination and Degradation of Bcr-Abl Protein

Daniel Alvira, Ruth Naughton, Lavinia Bhatt, Sara Tedesco, William D. Landry, and Thomas G. Cotter
J. Biol. Chem., Sep 2011; 286: 32313 - 32323.

Proteases

Doxycycline, a matrix metalloprotease inhibitor, reduces vascular remodeling and damage after cerebral ischemia in stroke-prone spontaneously hypertensive rats

Paulo W. Pires, Curt T. Rogers, Jonathon L. McClain, Hannah S. Garver, Gregory D. Fink, and Anne M. Dorrance
Am J Physiol Heart Circ Physiol, Jul 2011; 301: H87 - H97.

Yapsin 1 immunoreactivity in α -cells of human pancreatic islets: implications for the processing of human proglucagon by mammalian aspartic proteases

Niamh X Cawley, Guida Portela-Gomes, Hong Lou, and Y Peng Loh
J. Endocrinol., Aug 2011; 210: 181 - 187.

Novel Positron Emission Tomography Tracer Distinguishes Normal from Cancerous Cells

Muhammad Saeed, David Sheff, and Amnon Kohen
J. Biol. Chem., Sep 2011; 286: 33872 - 33878.

Neural circuit architecture defects in a *Drosophila* model of Fragile X syndrome are alleviated by minocycline treatment and genetic removal of matrix metalloproteinase

Saul S. Siller and Kendal Broadie
Dis. Model. Mech., Sep 2011; 4: 673 - 685.

Receptors

Chemerin, a Novel Peroxisome Proliferator-activated Receptor γ (PPAR γ) Target Gene That Promotes Mesenchymal Stem Cell Adipogenesis

Shanmugam Muruganandan, Sebastian D. Parlee, Jillian L. Rourke, Matthew C. Ernst, Kerry B. Goralski, and Christopher J. Sinal
J. Biol. Chem., Jul 2011; 286: 23982 - 23995.

Occupation of Either Site for the Neurosteroid Allopregnanolone Potentiates the Opening of the GABA_A Receptor Induced from Either Transmitter Binding Site

John Bracamontes, Megan McCollum, Caroline Esch, Ping Li, Jason Ann, Joe Henry Steinbach, and Gustav Akk
Mol. Pharmacol., Jul 2011; 80: 79 - 86.

Thioredoxin-Interacting Protein Mediates TRX1 Translocation to the Plasma Membrane in Response to Tumor Necrosis Factor- α : A Key Mechanism for Vascular Endothelial Growth Factor Receptor-2 Transactivation by Reactive Oxygen Species

Cameron World, Oded N. Spindel, and Bradford C. Berk
Arterioscler Thromb Vasc Biol, Aug 2011; 31: 1890 - 1897.

Nuclear Receptors TR2 and TR4 Recruit Multiple Epigenetic Transcriptional Corepressors That Associate Specifically with the Embryonic β -Type Globin Promoters in Differentiated Adult Erythroid Cells

Shuaiying Cui, Katarzyna E. Kolodziej, Naoshi Obara, Alexandra Amaral-Psarris, Jeroen Demmers, Lihong Shi, James Douglas Engel, Frank Grosveld, John Strouboulis, and Osamu Tanabe
Mol. Cell. Biol., Aug 2011; 31: 3298 - 3311.

NMDA Receptor Signaling in Oligodendrocyte Progenitors Is Not Required for Oligodendrogenesis and Myelination

Lindsay M. De Biase, Shin H. Kang, Emily G. Baxi, Masahiro Fukaya, Michele L. Pucak, Masayoshi Mishina, Peter A. Calabresi, and Dwight E. Bergles
J. Neurosci., Aug 2011; 31: 12650 - 12662.

G Protein Binding Sites on Calnuc (Nucleobindin 1) and NUCB2 (Nucleobindin 2) Define a New Class of G α_i -regulatory Motifs

Mikel Garcia-Marcos, Patrick S. Kietsunthorn, Honghui Wang, Pradipta Ghosh, and Marilyn G. Farquhar
J. Biol. Chem., Aug 2011; 286: 28138 - 28149.

Engineered Bivalent Ligands to Bias ErbB Receptor-mediated Signaling and Phenotypes

Steven M. Jay, Elma Kurtagic, Luis M. Alvarez, Seymour de Picciotto, Edgar Sanchez, Jessica F. Hawkins, Robin N. Prince, Yadir Guerrero, Carolyn L. Treasure, Richard T. Lee, and Linda G. Griffith
J. Biol. Chem., Aug 2011; 286: 27729 - 27740.

Characterization of a Sphingosine 1-Phosphate Receptor Antagonist Prodrug

Perry C. Kennedy, Ran Zhu, Tao Huang, Jose L. Tomsig, Thomas P. Mathews, Marion David, Olivier Peyruchaud, Timothy L. Macdonald, and Kevin R. Lynch
J. Pharmacol. Exp. Ther., Sep 2011; 338: 879 - 889.

Cytokine-Induced Alterations of $\alpha 7$ Nicotinic Receptor in Colonic CD4 T Cells Mediate Dichotomous Response to Nicotine in Murine Models of Th1/Th17- versus Th2-Mediated Colitis

Valentin Galitovskiy, Jing Qian, Alexander I. Chernyavsky, Steve Marchenko, Vivian Gindi, Robert A. Edwards, and Sergei A. Grando
J. Immunol., Sep 2011; 187: 2677 – 2687.

Role of Key Aromatic Residues in the Ligand-binding Domain of $\alpha 7$ Nicotinic Receptors in the Agonist Action of β -Amyloid

Mei Tong, Komal Arora, Michael M. White, and Robert A. Nichols
J. Biol. Chem., Sep 2011; 286: 34373 - 34381.

The Cytosolic Chaperone Hsc70 Promotes Traffic to the Cell Surface of Intracellular Retained Melanocortin-4 Receptor Mutants

Eirini Meimaridou, Sakina B. Gooljar, Nalini Ramnarace, Lucia Anthonypillai, Adrian J. L. Clark, and J. Paul Chapple
Mol. Endocrinol., Sep 2011; 25: 1650 - 1660.

A selective estrogen receptor α agonist ameliorates hepatic steatosis in the male aromatase knockout mouse

Jenny DY Chow, Margaret E E Jones, Katja Prella, Evan R Simpson, and Wah Chin Boon
J. Endocrinol., Sep 2011; 210: 323 - 334.

RNAi

The Tpr protein regulates export of mRNAs with retained introns that traffic through the Nxf1 pathway

John H. Coyle, Yeou-Cherng Bor, David Rekosh, and Marie-Louise Hammarskjold
RNA, Jul 2011; 17: 1344 - 1356.

Chloride Channel (Clc)-5 Is Necessary for Exocytic Trafficking of Na⁺/H⁺ Exchanger 3 (NHE3)

Zhihong Lin, Shi Jin, Xiaohong Duan, Tong Wang, Sabrina Martini, Phuson Hulamm, Boyoung Cha, Ann Hubbard, Mark Donowitz, and Sandra E. Guggino
J. Biol. Chem., Jul 2011; 286: 22833 - 22845.

The histone methyltransferase Set7/9 promotes myoblast differentiation and myofibril assembly

Yazhong Tao, Ronald L. Nepl, Zhan-Peng Huang, Jianfu Chen, Ru-Hang Tang, Ru Cao, Yi Zhang, Suk-Won Jin, and Da-Zhi Wang
J. Cell Biol., Aug 2011; 194: 551 - 565.

Dual Function of Protein Kinase C (PKC) in 12-O-Tetradecanoylphorbol-13-acetate (TPA)-induced Manganese Superoxide Dismutase (MnSOD) Expression: ACTIVATION OF CREB AND FOXO3a BY PKC- α PHOSPHORYLATION AND BY PKC-MEDIATED INACTIVATION OF Akt, RESPECTIVELY

Youn Wook Chung, Ha Kun Kim, IckYoung Kim, Moon B. Yim, and P. Boon Chock
J. Biol. Chem., Aug 2011; 286: 29681 - 29690.

Development of a Highly Sensitive Cytotoxicity Assay System for CYP3A4-Mediated Metabolic Activation

Hiroko Hosomi, Tatsuki Fukami, Atsushi Iwamura, Miki Nakajima, and Tsuyoshi Yokoi
Drug Metab. Dispos., Aug 2011; 39: 1388 - 1395.

APOBEC3G Complexes Decrease Human Immunodeficiency Virus Type 1 Production

Kenneth L. Martin, Megan Johnson, and Richard T. D'Aquila
J. Virol., Sep 2011; 85: 9314 - 9326.

The Protein Kinase IKK ϵ Is a Potential Target for the Treatment of Inflammatory Hyperalgesia

Christine V. Möser, Katharina Kynast, Katharina Baatz, Otto Quintus Russe, Nerea Ferreirós, Heike Costiuk, Ruirui Lu, Achim Schmidtke, Irmgard Tegeder, Gerd Geisslinger, and Ellen Niederberger
J. Immunol., Sep 2011; 187: 2617 - 2625.

Inhibition of Protein-tyrosine Phosphatase 1B (PTP1B) Mediates Ubiquitination and Degradation of Bcr-Abl Protein

Daniel Alvira, Ruth Naughton, Lavinia Bhatt, Sara Tedesco, William D. Landry, and Thomas G. Cotter
J. Biol. Chem., Sep 2011; 286: 32313 - 32323.

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., Sep 2011; 20: 3356 - 3365.

ROS – Reactive Oxygen Species

E4F1 deficiency results in oxidative stress-mediated cell death of leukemic cells

Elodie Hatchi, Genevieve Rodier, Matthieu Lacroix, Julie Caramel, Olivier Kirsh, Chantal Jacquet, Emilie Schrepfer, Sylviane Lagarrigue, Laetitia Karine Linares, Gwendaline Lledo, Sylvie Tondeur, Pierre Dubus, Claude Sardet, and Laurent Le Cam
J. Exp. Med., Jul 2011; 208: 1403 - 1417.

GSH monoethyl ester rescues mitochondrial defects in cystic fibrosis models

Mairead Kelly-Aubert, Stéphanie Trudel, Janine Fritsch, Thao Nguyen-Khoa, Maryvonne Baudouin-Legros, Sandra Moriceau, Ludovic Jeanson, Fatima Djouadi, Corine Matar, Marc Conti, Mario Ollero, Franck Brouillard, and Aleksander Edelman
Hum. Mol. Genet., Jul 2011; 20: 2745 - 2759.

Regulation of the expression of the avian uncoupling protein 3 by isoproterenol and fatty acids in chick myoblasts: possible involvement of AMPK and PPAR α ?

Romain Joubert, Sonia Métayer-Coustard, Sabine Crochet, Estelle Cailleau-Audouin, Joëlle Dupont, Michel Jacques Duclos, Sophie Tesseraud, and Anne Collin
Am J Physiol Regulatory Integrative Comp Physiol, Jul 2011; 301: R201 - R208.

Thioredoxin-Interacting Protein Mediates TRX1 Translocation to the Plasma Membrane in Response to Tumor Necrosis Factor- α : A Key Mechanism for Vascular Endothelial Growth Factor Receptor-2 Transactivation by Reactive Oxygen Species

Cameron World, Oded N. Spindel, and Bradford C. Berk
Arterioscler Thromb Vasc Biol, Aug 2011; 31: 1890 - 1897.

Dual Function of Protein Kinase C (PKC) in 12-O-Tetradecanoylphorbol-13-acetate (TPA)-induced Manganese Superoxide Dismutase (MnSOD) Expression: ACTIVATION OF CREB AND FOXO3a BY PKC- α PHOSPHORYLATION AND BY PKC-MEDIATED INACTIVATION OF Akt, RESPECTIVELY

Youn Wook Chung, Ha Kun Kim, Ick Young Kim, Moon B. Yim, and P. Boon Chock
J. Biol. Chem., Aug 2011; 286: 29681 - 29690.

Inhibition of Protein-tyrosine Phosphatase 1B (PTP1B) Mediates Ubiquitination and Degradation of Bcr-Abl Protein

Daniel Alvira, Ruth Naughton, Lavinia Bhatt, Sara Tedesco, William D. Landry, and Thomas G. Cotter
J. Biol. Chem., Sep 2011; 286: 32313 - 32323.

Transcription Factors

Thioredoxin-Interacting Protein Mediates TRX1 Translocation to the Plasma Membrane in Response to Tumor Necrosis Factor- α : A Key Mechanism for Vascular Endothelial Growth Factor Receptor-2 Transactivation by Reactive Oxygen Species

Cameron World, Oded N. Spindel, and Bradford C. Berk
Arterioscler Thromb Vasc Biol, Aug 2011; 31: 1890 - 1897.

Simvastatin Inhibits IFN Regulatory Factor 4 Expression and Th17 Cell Differentiation in CD4⁺T Cells Derived from Patients with Multiple Sclerosis

Xin Zhang, Yazhong Tao, Luigi Troiani, and Silva Markovic-Plese
J. Immunol., Sep 2011; 187: 3431 - 3437.

Ubiquitination and Sumoylation

SUMO conjugation of STAT1 protects cells from hyperresponsiveness to IFN γ

Andreas Begitt, Mathias Droscher, Klaus-Peter Knobeloch, and Uwe Vinkemeier
Blood, Jul 2011; 118: 1002 - 1007.

The E3 ubiquitin ligase specificity subunit ASB2 α targets filamins for proteasomal degradation by interacting with the filamin actin-binding domain

Ziba Razinia, Massimiliano Baldassarre, Mohamed Bouaouina, Isabelle Lamsoul, Pierre G. Lutz, and David A. Calderwood
J. Cell Sci., Aug 2011; 124: 2631 - 2641.

Inhibition of Protein-tyrosine Phosphatase 1B (PTP1B) Mediates Ubiquitination and Degradation of Bcr-Abl Protein

Daniel Alvira, Ruth Naughton, Lavinia Bhatt, Sara Tedesco, William D. Landry, and Thomas G. Cotter
J. Biol. Chem., Sep 2011; 286: 32313 - 32323.

Agonist-directed Interactions with Specific β -Arrestins Determine μ -Opioid Receptor Trafficking, Ubiquitination, and Dephosphorylation

Chad E. Groer, Cullen L. Schmid, Alex M. Jaeger, and Laura M. Bohn
J. Biol. Chem., Sep 2011; 286: 31731 - 31741.

Cleavage of the Adaptor Protein TRIF by Enterovirus 71 3C Inhibits Antiviral Responses Mediated by Toll-Like Receptor 3

Xiaobo Lei, Zhenmin Sun, Xinlei Liu, Qi Jin, Bin He, and Jianwei Wang
J. Virol., Sep 2011; 85: 8811 - 8818.

WNT Protein-independent Constitutive Nuclear Localization of β -Catenin Protein and Its Low Degradation Rate in Thalamic Neurons

Katarzyna Misztal, Marta B. Wisniewska, Mateusz Ambrozkiwicz, Andrzej Nagalski, and Jacek Kuznicki
J. Biol. Chem., Sep 2011; 286: 31781 - 31788.

Mono-ubiquitination Drives Nuclear Export of the Human DCN1-like Protein hDCNL1

Kenneth Wu, Hua Yan, Lei Fang, Xinjiang Wang, Cathie Pflieger, Xuejun Jiang, Lan Huang, and Zhen-Qiang Pan
J. Biol. Chem., Sep 2011; 286: 34060 - 34070.

MARCH ubiquitin ligases alter the itinerary of clathrin-independent cargo from recycling to degradation

Craig A. Eyster, Nelson B. Cole, Shariska Petersen, Kasinath Viswanathan, Klaus Früh, and Julie G. Donaldson
Mol. Biol. Cell, Sep 2011; 22: 3218 - 3230.

Virology/Microbiology

Influence of PB2 host-range determinants on the intranuclear mobility of the influenza A virus polymerase

Ágnes Foeglein, Eva M. Loucaides, Manuela Mura, Helen M. Wise, Wendy S. Barclay, and Paul Digard
J. Gen. Virol., Jul 2011; 92: 1650 - 1661.

TREM-1 Amplifies Corneal Inflammation after *Pseudomonas aeruginosa* Infection by Modulating Toll-Like Receptor Signaling and Th1/Th2-Type Immune Responses

Minhao Wu, Anping Peng, Mingxia Sun, Qiuchan Deng, Linda D. Hazlett, Jin Yuan, Xialin Liu, Qianying Gao, Lianqiang Feng, Junfang He, Ping Zhang, and Xi Huang
Infect. Immun., Jul 2011; 79: 2709 - 2716.

Placental Expression of the Heme Transporter, Feline Leukemia Virus Subgroup C Receptor, Is related to Maternal Iron Status in Pregnant Adolescents

Lindsay M. Jaacks, Melissa F. Young, Bridget V. Essley, Thomas J. McNanley, Elizabeth M. Cooper, Eva K. Pressman, Allison W. McIntyre, Mark S. Orlando, Janis L. Abkowitz, Ronnie Guillet, and Kimberly O. O'Brien
J. Nutr., Jul 2011; 141: 1267 - 1272.

The PI3K/Akt pathway is involved in early infection of some exogenous avian leukosis viruses

Shao-zhen Feng, Wei-sheng Cao, and Ming Liao
J. Gen. Virol., Jul 2011; 92: 1688 - 1697.

Human Immunodeficiency Virus Type 1 Modified To Package Simian Immunodeficiency Virus Vpx Efficiently Infects Macrophages and Dendritic Cells

Nicole Sunseri, Meagan O'Brien, Nina Bhardwaj, and Nathaniel R. Landau
J. Virol., Jul 2011; 85: 6263 - 6274.

Sprouty 2 Binds ESCRT-II Factor Eap20 and Facilitates HIV-1 Gag Release

G. N. Medina, L. S. Ehrlich, M. H. Chen, M. B. Khan, M. D. Powell, and C. A. Carter
J. Virol., Jul 2011; 85: 7353 - 7362.

Protein Kinase C and NF- κ B-Dependent CD4 Downregulation in Macrophages Induced by T Cell-Derived Soluble Factors: Consequences for HIV-1 Infection

Rui André Saraiva Raposo, David C. Trudgian, Benjamin Thomas, Bonnie van Wilgenburg, Sally A. Cowley, and William James
J. Immunol., Jul 2011; 187: 748 - 759.

Absence of XMRV Retrovirus and Other Murine Leukemia Virus-Related Viruses in Patients with Chronic Fatigue Syndrome

Clifford H. Shin, Lucinda Bateman, Robert Schlaberg, Ashley M. Bunker, Christopher J. Leonard, Ronald W. Hughen, Alan R. Light, Kathleen C. Light, and Ila R. Singh
J. Virol., Jul 2011; 85: 7195 - 7202.

Hepatitis C Virus Infection Causes Cell Cycle Arrest at the Level of Initiation of Mitosis

Rathi P. Kannan, Lucinda L. Hensley, Lauren E. Evers, Stanley M. Lemon, and David R. McGivern
J. Virol., Aug 2011; 85: 7989 - 8001.

Analysis of Rab GTPase-Activating Proteins Indicates that Rab1a/b and Rab43 Are Important for Herpes Simplex Virus 1 Secondary Envelopment

Helen L. Zenner, Shin-ichiro Yoshimura, Francis A. Barr, and Colin M. Crump
J. Virol., Aug 2011; 85: 8012 - 8021.

Classical Swine Fever Virus N^{pro} Limits Type I Interferon Induction in Plasmacytoid Dendritic Cells by Interacting with Interferon Regulatory Factor 7

Ana R. Fiebach, Laurence Guzylack-Piriou, Sylvie Python, Artur Summerfield, and Nicolas Ruggli
J. Virol., Aug 2011; 85: 8002 - 8011.

Host Response to Polyomavirus Infection Is Modulated by RNA Adenosine Deaminase ADAR1 but Not by ADAR2

Cyril X. George and Charles E. Samuel
J. Virol., Aug 2011; 85: 8338 - 8347.

The Cytoplasmic Domain of Marburg Virus GP Modulates Early Steps of Viral Infection

Eva Mittler, Larissa Kolesnikova, Bettina Hartlieb, Robert Davey, and Stephan Becker
J. Virol., Aug 2011; 85: 8188 - 8196.

The Herpes Simplex Virus 1 UL17 Protein Is the Second Constituent of the Capsid Vertex-Specific Component Required for DNA Packaging and Retention

Katerina Toropova, Jamie B. Huffman, Fred L. Homa, and James F. Conway
J. Virol., Aug 2011; 85: 7513 - 7522.

Herpes Simplex Virus Immediate-Early Protein ICP0 Is Targeted by SIAH-1 for Proteasomal Degradation

Claus-Henning Nagel, Nina Albrecht, Kristijana Milovic-Holm, Lakshmikanth Mariyanna, Britta Keyser, Bettina Abel, Britta Weseloh, Thomas G. Hofmann, Martha M. Eibl, and Joachim Hauber
J. Virol., Aug 2011; 85: 7644 - 7657.

Hepatitis C Virus Infection Causes Cell Cycle Arrest at the Level of Initiation of Mitosis

Rathi P. Kannan, Lucinda L. Hensley, Lauren E. Evers, Stanley M. Lemon, and David R. McGivern
J. Virol., Aug 2011; 85: 7989 - 8001.

Hepatitis C Virus (HCV) Envelope Glycoproteins E1 and E2 Contain Reduced Cysteine Residues Essential for Virus Entry

Johanna Fraser, Irene Boo, Pantelis Pountourios, and Heidi E. Drummer
J. Biol. Chem., Sep 2011; 286: 31984 - 31992.

Classical Swine Fever Virus N^{pro} Limits Type I Interferon Induction in Plasmacytoid Dendritic Cells by Interacting with Interferon Regulatory Factor 7

Ana R. Fiebach, Laurence Guzylack-Piriou, Sylvie Python, Artur Summerfield, and Nicolas Ruggli
J. Virol., Aug 2011; 85: 8002 - 8011.

Daedalus: a robust, turnkey platform for rapid production of decigram quantities of active recombinant proteins in human cell lines using novel lentiviral vectors

Ashok D. Bandaranayake, Colin Correnti, Byoung Y. Ryu, Michelle Brault, Roland K. Strong, and David J. Rawlings
Nucleic Acids Res., Sep 2011; 10.1093/nar/gkr706.

APOBEC3G Complexes Decrease Human Immunodeficiency Virus Type 1 Production

Kenneth L. Martin, Megan Johnson, and Richard T. D'Aquila
J. Virol., Sep 2011; 85: 9314 - 9326.

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., Sep 2011; 39: 7775 - 7790

APPLICATIONS

In-Cell Western™ Assays

Endorepellin, the Angiostatic Module of Perlecan, Interacts with Both the $\alpha 2\beta 1$ Integrin and Vascular Endothelial Growth Factor Receptor 2 (VEGFR2): A DUAL RECEPTOR ANTAGONISM

Atul Goyal, Nutan Pal, Matthew Concannon, Matthew Paul, Mike Doran, Chiara Poluzzi, Kiyotoshi Sekiguchi, John M. Whitelock, Thomas Neill, and Renato V. Iozzo
J. Biol. Chem., Jul 2011; 286: 25947 - 25962.

CD82 gene suppression in endometrial stromal cells leads to increase of the cell invasiveness in the endometriotic milieu

Ming-Qing Li, Xiao-Fan Hou, Shi-Jian Lv, Yu-Han Meng, Xiao-Qiu Wang, Chuan-Ling Tang, and Da-Jin Li
J. Mol. Endocrinol., Aug 2011; 47: 195 - 208.

Erythropoietin Increases Expression and Function of Transient Receptor Potential Canonical 5 Channels

Ying Liu, Yunfei Xu, Florian Thilo, Ulla G. Friis, Boye L. Jensen, Alexandra Scholze, Junhua Zheng, and Martin Tepel
Hypertension, Aug 2011; 58: 317 - 324.

Cytokine-Induced Alterations of $\alpha 7$ Nicotinic Receptor in Colonic CD4 T Cells Mediate Dichotomous Response to Nicotine in Murine Models of Th1/Th17- versus Th2-Mediated Colitis

Valentin Galitovskiy, Jing Qian, Alexander I. Chernyavsky, Steve Marchenko, Vivian Gindi, Robert A. Edwards, and Sergei A. Grando
J. Immunol., Sep 2011; 187: 2677 - 2687.

The Cytosolic Chaperone Hsc70 Promotes Traffic to the Cell Surface of Intracellular Retained Melanocortin-4 Receptor Mutants

Eirini Meimaridou, Sakina B. Gooljar, Nalini Ramnarace, Lucia Anthonypillai, Adrian J. L. Clark, and J. Paul Chapple
Mol. Endocrinol., Sep 2011; 25: 1650 - 1660.

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® Accessory

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, Andréas Bikfalvi, and Hervé 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 (8)

BIOCHEMISTRY:

Unveiling the structural basis for translational ambiguity tolerance in a human fungal pathogen

Rita Rocha, Pedro José Barbosa Pereira, Manuel A. S. Santos, and Sandra Macedo-Ribeiro
PNAS, Aug 2011; 108: 14091 - 14096.

Processive phosphorylation of ERK MAP kinase in mammalian cells

Kazuhiro Aoki, Masashi Yamada, Katsuyuki Kunida, Shuhei Yasuda, and Michiyuki Matsuda
PNAS, Aug 2011; 108: 12675 - 12680.

CELL BIOLOGY:

From the Cover: Feature Article: A haploid genetic screen identifies the major facilitator domain containing 2A (MFSD2A) transporter as a key mediator in the response to tunicamycin

Jan H. Reiling, Clary B. Clish, Jan E. Carette, Malini Varadarajan, Thijn R. Brummelkamp, and David M. Sabatini
PNAS, Jul 2011; 108: 11756 - 11765.

ENGINEERING:

Blast-induced phenotypic switching in cerebral vasospasm

Patrick W. Alford, Borna E. Dabiri, Josue A. Goss, Matthew A. Hemphill, Mark D. Brigham, and Kevin Kit Parker
PNAS, Aug 2011; 108: 12705 - 12710.

GENETICS:

Neuronal circuitry regulates the response of *Caenorhabditis elegans* to misfolded proteins

Veena Prahlad and Richard I. Morimoto
PNAS, Aug 2011; 108: 14204 - 14209.

NEUROSCIENCE:

CNS expression of glucocerebrosidase corrects α -synuclein pathology and memory in a mouse model of Gaucher-related synucleinopathy

S. Pablo Sardi, Jennifer Clarke, Cathrine Kinnecom, Thomas J. Tamsett, Lingyun Li, Lisa M. Stanek, Marco A. Passini, Gregory A. Grabowski, Michael G. Schlossmacher, Richard L. Sidman, Seng H. Cheng, and Lamya S. Shihabuddin
PNAS, Jul 2011; 108: 12101 - 12106.

S-Nitrosylation activates Cdk5 and contributes to synaptic spine loss induced by β -amyloid peptide

Jing Qu, Tomohiro Nakamura, Gang Cao, Emily A. Holland, Scott R. McKercher, and Stuart A. Lipton
PNAS, Aug 2011; 108: 14330 - 14335.

The neuronal K-Cl cotransporter KCC2 influences postsynaptic AMPA receptor content and lateral diffusion in dendritic spines

Grégory Gouvain, Ingrid Chamma, Quentin Chevy, Carolina Cabezas, Theano Irinopoulou, Natalia Bodrug, Michèle Carnaud, Sabine Lévi, and Jean Christophe Poncer
PNAS, Sep 2011; 108: 15474 - 15479.

©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 I 1011