

Recent Odyssey® Publications

(Selected from 285 publications in June 09 – August 09 HighWire Search)

Fall 2009

Volume 5

www.highwire.org currently references more than 2,200 peer-reviewed publications using the LI-COR® Odyssey® Infrared Imaging System. The publications listed in this issue highlight the wide range of research areas and applications in which the Odyssey is used.

TABLE OF CONTENTS

	Page
Apoptosis2
Angiogenesis2
Autophagy2
Cancer3
Cardiovascular3-4
Cell Adhesion, Migration and Invasion4
Cell Signaling4-5
Cell Cycle5-6
Hypoxia6
Ion Channels6
Metabolism6-7
Matrix Metaloproteases7
Muscle7
Neuroscience7-8
Nephrology8
Phosphatases8
Receptors8-9
RNAi9
Stem Cells9
Transcription Factors10
Ubiquitination and Sumoylation10
Virology11
Applications11-12
<i>In-Cell Western™</i>	
<i>Quantitative Western (Molecules/Cell)</i>	
<i>MousePod</i>	
<i>Quantitative Reverse Phase Protein Arrays</i>	
<i>DNA Methylation Assay</i>	

PUBLICATIONS REFERENCING ODYSSEY INFRARED IMAGING DATA

RESEARCH AREAS

APOPTOSIS

Bufalin Induces Reactive Oxygen Species Dependent Bax Translocation and Apoptosis in ASTC-a-1 Cells

Lei Sun, Tongsheng Chen, Xiaoping Wang, Yun Chen, and Xunbin Wei
Evid. Based Complement. Altern. Med., Jul 2009; 10.1093/ecam/nep082.

CMTM3, Located at the Critical Tumor Suppressor Locus 16q22.1, Is Silenced by CpG Methylation in Carcinomas and Inhibits Tumor Cell Growth through Inducing Apoptosis

Yu Wang, Jisheng Li, Yan Cui, Ting Li, Ka Man Ng, Hua Geng, Henan Li, Xing-sheng Shu, Hongyu Li, Wei Liu, Bing Luo, Qian Zhang, Tony Shu Kam Mok, Wei Zheng, Xiaoyan Qiu, Gopesh Srivastava, Jun Yu, Joseph J.Y. Sung, Anthony T.C. Chan, Dalong Ma, Qian Tao, and Wenling Han
Cancer Res., Jun 2009; 69: 5194 - 5201

PUMA Promotes Bax Translocation by Both Directly Interacting with Bax and by Competitive Binding to Bcl-XL during UV-induced Apoptosis

Yingjie Zhang, Da Xing, and Lei Liu
Mol. Biol. Cell, Jul 2009; 20: 3077 - 3087.

Rab7 Activation by Growth Factor Withdrawal Contributes to the Induction of Apoptosis

Kimberly Romero Rosales, Eigen R. Peralta, Garret G. Guenther, Susan Y. Wong, and Aimee L. Edinger
Mol. Biol. Cell, Jun 2009; 20: 2831 - 2840.

Real-Time Detection of Caspase-3-Like Protease Activation *in Vivo* Using Fluorescence Resonance Energy Transfer during Plant Programmed Cell Death Induced by Ultraviolet C Overexposure

Lingrui Zhang, Qixian Xu, Da Xing, Caiji Gao, and Hongwu Xiong
Plant Physiology, Aug 2009; 150: 1773 - 1783.

Serine 58 of 14-3-3 Is a Molecular Switch Regulating ASK1 and Oxidant Stress-Induced Cell Death

Jibin Zhou, Zhili Shao, Risto Kerkela, Hidenori Ichijo, Anthony J. Muslin, Celia Pombo, and Thomas Force
Mol. Cell. Biol., Aug 2009; 29: 4167 - 4176.

ANGIOGENESIS

Phosphorylation of endothelial nitric oxide synthase by atypical PKC contributes to angiotensin-1-dependent inhibition of VEGF-induced endothelial permeability *in vitro*

Malika Oubaha and Jean-Philippe Gratton
Blood, Jun 2009; 10.1182/blood-2008-12-196584.

Pathogenic role of angiotensin II and oxidized LDL in obstructive sleep apnoea

T. Kizawa, Y. Nakamura, S. Takahashi, S. Sakurai, K. Yamauchi, and H. Inoue
Eur. Respir. J., Jul 2009; 10.1183/09031936.00009709.

AUTOPHAGY

Degradation of regulator of calcineurin 1 (RCAN1) is mediated by both chaperone-mediated autophagy and ubiquitin proteasome pathways

Heng Liu, Pin Wang, Weihong Song, and Xiulian Sun
FASEB J, Jun 2009; 10.1096/fj.09-134296.

Clathrin Functions in the Absence of the Terminal Domain Binding Site for Adaptor-associated Clathrin-Box Motifs

John R. Collette, Richard J. Chi, Douglas R. Boettner, Isabel M. Fernandez-Golbano, Rachael Plemel, Alex J. Merz, Maria Isabel Geli, Linton M. Traub, and Sandra K. Lemmon
Mol. Biol. Cell, Jul 2009; 20: 3401 - 3413.

CANCER

EGFRvIII and c-Met pathway inhibitors synergize against PTEN-null/EGFRvIII+ glioblastoma xenografts

Bachchu Lal, C. Rory Goodwin, Yingying Sang, Catherine A. Foss, Kathrine Cornet, Sameena Muzamil, Martin G. Pomper, Jin Kim, and John Laterra
Mol. Cancer Ther., Jul 2009; 8: 1751 - 1760.

Activation of Rap1 Promotes Prostate Cancer Metastasis

Candice L. Bailey, Patrick Kelly, and Patrick J. Casey
Cancer Res., Jun 2009; 69: 4962 - 4968.

Novel mechanism of IGF-binding protein-3 action on prostate cancer cells: inhibition of proliferation, adhesion, and motility

Petra Massoner, Daniela Colleselli, Andrea Matscheski, Haymo Pircher, Stephan Geley, Pidder Jansen Dürr, and Helmut Klocker
Endocr. Relat. Cancer, Sep 2009; 16: 795 - 808.

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.

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., Jul 2009; 69: 5575 - 5583.

Targeting Focal Adhesion Kinase with Dominant-Negative FRNK or Hsp90 Inhibitor 17-DMAG Suppresses Tumor Growth and Metastasis of SiHa Cervical Xenografts

Joerg Schwock, Neesha Dhani, Mary Ping-Jiang Cao, Jinzi Zheng, Richard Clarkson, Nikolina Radulovich, Roya Navab, Lars-Christian Horn, and David W. Hedley
Cancer Res., Jun 2009; 69: 4750 - 4759.

Sorafenib and Vorinostat Kill Colon Cancer Cells by CD95-Dependent and -Independent Mechanisms

Teneille Walker, Clint Mitchell, Margaret A. Park, Adly Yacoub, Martin Graf, Mohamed Rahmani, Peter J. Houghton, Christina Voelkel-Johnson, Steven Grant, and Paul Dent
Mol. Pharmacol., Aug 2009; 76: 342 - 355.

The tumor suppressor functions of p27^{kip1} include the control of mesenchymal/amoeboid transition

Stefania Berton, Barbara Belletti, Katarina Wolf, Vincenzo Canzonieri, Francesca Lovat, Andrea Vecchione, Alfonso Colombatti, Peter Friedl, and Gustavo Baldassarre
Mol. Cell. Biol., Jul 2009; 10.1128/MCB.00144-09.

Novel Mechanism of Lapatinib Resistance in HER2-Positive Breast Tumor Cells: Activation of AXL

Li Liu, James Greger, Hong Shi, Yuan Liu, Joel Greshock, Roland Annan, Wendy Halsey, Ganesh M. Sathe, Anne-Marie Martin, and Tona M. Gilmer
Cancer Res., Aug 2009; 10.1158/0008-5472.CAN-08-4490.

CARDIOVASCULAR

Phenylephrine and sustained acidosis activate the neonatal rat cardiomyocyte Na⁺/H⁺ exchanger through phosphorylation of amino acids Ser⁷⁷⁰ and Ser⁷⁷¹

Ersilia Coccaro, Pratap Karki, Cicerone Cojocar, and Larry Fliegel
Am J Physiol Heart Circ Physiol, Aug 2009; 297: H846 - H858.

TRPC1 and STIM1 mediate capacitative Ca²⁺ entry in mouse pulmonary arterial smooth muscle cells

Lih Chyuan Ng, Mary D. McCormack, Judith A. Airey, Cherie A. Singer, Phillip S. Keller, Xiao-Ming Shen, and Joseph R. Hume
J. Physiol., Jun 2009; 587: 2429 - 2442.

Hindlimb Unweighting Induces Changes in the p38^{MAPK} Contractile Pathway of the Rat Abdominal Aorta

Scott Matthew Summers, Yuichiro Hayashi, Steven Vu Nguyen, Thu Minh Nguyen, and Ralph Earl Purdy
J Appl Physiol, July 2009; 107: 121 - 127.

CELL ADHESION, MIGRATION AND INVASION

Splicing factors PTBP1 and PTBP2 promote proliferation and migration of glioma cell lines

Hannah C. Cheung, Tao Hai, Wen Zhu, Keith A. Baggerly, Spiridon Tsavachidis, Ralf Krahe, and Gilbert J. Cote
Brain, Aug 2009; 132: 2277 - 2288.

The Pyk2 FERM domain as a target to inhibit glioma migration

Joseph C. Loftus, Zhongbo Yang, Nhan L. Tran, Jean Kloss, Carole Viso, Michael E. Berens, and Christopher A. Lipinski
Mol. Cancer Ther., Jun 2009; 8: 1505 - 1514.

Phosphorylation of RACK1 on Tyrosine 52 by c-Abl Is Required for Insulin-like Growth Factor I-mediated Regulation of Focal Adhesion Kinase

Patrick A. Kiely, George S. Baillie, Robert Barrett, Deirdre A. Buckley, David R. Adams, Miles D. Houslay, and Rosemary O'Connor
J. Biol. Chem., Jul 2009; 284: 20263 - 20274.

Up-regulation of miR-21 by HER2/*neu* Signaling Promotes Cell Invasion

Tzu-Hsuan Huang, Fangting Wu, Gabriel B. Loeb, Ruby Hsu, Amy Heidersbach, Allison Brincat, Dai Horiuchi, Robert J. Lebbink, Yin-Yuan Mo, Andrei Goga, and Michael T. McManus
J. Biol. Chem., Jul 2009; 284: 18515 - 18524.

Breast Cancer Migration and Invasion Depend on Proteasome Degradation of Regulator of G-Protein Signaling 4

Yan Xie, Dennis W. Wolff, Taotao Wei, Bo Wang, Caishu Deng, Joseph K. Kirui, Haihong Jiang, Jianbing Qin, Peter W. Abel, and Yaping Tu
Cancer Res., Jul 2009; 69: 5743 - 5751.

Combined Inhibition of MEK and Mammalian Target of Rapamycin Abolishes Phosphorylation of Cyclin-Dependent Kinase 4 in Glioblastoma Cell Lines and Prevents Their Proliferation

Sabine Paternot and Pierre P. Roger
Cancer Res., Jun 2009; 69: 4577 - 4581.

Stretch-stimulated glucose uptake in skeletal muscle is mediated by reactive oxygen species and p38 MAP-kinase

Melissa A. Chambers, Jennifer S. Moylan, Jeffrey D. Smith, Laurie J. Goodyear, and Michael B. Reid
J. Physiol., Jul 2009; 587: 3363 - 3373.

Plectin Isoform-dependent Regulation of Keratin-Integrin $\alpha6\beta4$ Anchorage via Ca²⁺/Calmodulin

Julius Kostan, Martin Gregor, Gernot Walko, and Gerhard Wiche
J. Biol. Chem., Jul 2009; 284: 18525 - 18536.

CELL SIGNALING

***Campylobacter jejuni*-Induced Activation of Dendritic Cells Involves Cooperative Signaling through Toll-Like Receptor 4 (TLR4)-MyD88 and TLR4-TRIF Axes**

Vijay A. K. Rathinam, Daniel M. Appledorn, Kathleen A. Hoag, Andrea Amalfitano, and Linda S. Mansfield
Infect. Immun., June 2009; 77: 2499 - 2507.

Inhibition of mTOR Signaling in Parkinson's Disease Prevents L-DOPA-Induced Dyskinesia

Emanuela Santini, Myriam Heiman, Paul Greengard, Emmanuel Valjent, and Gilberto Fisone
Sci. Signal., Jul 2009; 2: ra36.

Sprouty2 Association with B-Raf Is Regulated by Phosphorylation and Kinase Conformation

Suzanne C. Brady, Mathew L. Coleman, June Munro, Stephan M. Feller, Nicolas A. Morrice, and Michael F. Olson
Cancer Res., Aug 2009; 10.1158/0008-5472.CAN-08-4447.

Intracellular Delivery of a Cell-Penetrating SOCS1 that Targets IFN- γ Signaling

Antonio DiGiandomenico, Lukasz S. Wylezinski, and Jacek Hawiger
Sci. Signal., Jul 2009; 2: ra37.

PI3K γ Adaptor Subunits Define Coupling to Degranulation and Cell Motility by Distinct PtdIns(3,4,5)P $_3$ Pools in Mast Cells

Thomas Bohnacker, Romina Marone, Emilie Collmann, Ronan Calvez, Emilio Hirsch, and Matthias P. Wymann
Sci. Signal., Jun 2009; 2: ra27.

Pivotal Advance: CEACAM1 is a negative coreceptor for the B cell receptor and promotes CD19-mediated adhesion of B cells in a PI3K-dependent manner

Elizabeth O. Lobo, Zhifang Zhang, and John E. Shively
J. Leukoc. Biol., Aug 2009; 86: 205 - 218.

Endosomal Endothelin-converting Enzyme-1: A REGULATOR OF β -ARRESTIN-DEPENDENT ERK SIGNALING

Graeme S. Cottrell, Benjamin E. Padilla, Silvia Amadesi, Daniel P. Poole, Jane E. Murphy, Markus Hardt, Dirk Roosterman, Martin Steinhoff, and Nigel W. Bunnett
J. Biol. Chem., Aug 2009; 284: 22411 - 22425.

Regulation of Jak2 Function by Phosphorylation of Tyr $_{317}$ and Tyr $_{637}$ during Cytokine Signaling

Scott A. Robertson, Rositsa I. Koleva, Lawrence S. Argetsinger, Christin Carter-Su, Jarrod A. Marto, Edward P. Feener, and Martin G. Myers, Jr.
Mol. Cell. Biol., Jun 2009; 29: 3367 - 3378.

Regulation of Platelet-derived Growth Factor Receptor Function by Integrin-associated Cell Surface Transglutaminase

Evgeny A. Zemskov, Elena Loukinova, Irina Mikhailenko, Richard A. Coleman, Dudley K. Strickland, and Alexey M. Belkin
J. Biol. Chem., Jun 2009; 284: 16693 - 16703.

MEK/ERK-Mediated Phosphorylation of Bim Is Required to Ensure Survival of T and B Lymphocytes during Mitogenic Stimulation

Lorraine A. O'Reilly, Elizabeth A. Kruse, Hamsa Puthalakath, Priscilla N. Kelly, Thomas Kaufmann, David C. S. Huang, and Andreas Strasser
J. Immunol., Jul 2009; 183: 261 - 269.

SIK1 Couples LKB1 to p53-Dependent Anoikis and Suppresses Metastasis

Hailing Cheng, Pixu Liu, Zhigang C. Wang, Lihua Zou, Stephanie Santiago, Victoria Garbitt, Ole V. Gjoerup, J. Dirk Iglehart, Alexander Miron, Andrea L. Richardson, William C. Hahn, and Jean J. Zhao
Sci. Signal., Jul 2009; 2: ra35.

O-GlcNAc modification of insulin receptor substrate-1 (IRS-1) occurs in close proximity to multiple SH2 domain binding motifs

Amanda L. Klein, Mary N. Berkaw, Maria G. Buse, and Lauren E. Ball
Mol. Cell. Proteomics, Aug 2009; 10.1074/mcp.M900207-MCP200.

CELL CYCLE

Dietary flavonoid fisetin induces a forced exit from mitosis by targeting the mitotic spindle checkpoint

Anna-Leena Salmela, Jeroen Pouwels, Asta Varis, Anu M. Kukkonen, Pauliina Toivonen, Pasi K. Halonen, Merja Perälä, Olli Kallioniemi, Gary J. Gorbisky, and Marko J. Kallio
Carcinogenesis, Jun 2009; 30: 1032 - 1040.

Effects of Nickel on Cyclin Expression, Cell Cycle Progression and Cell Proliferation in Human Pulmonary Cells

Jin Ding, Guoping He, Wenfeng Gong, Wen Wen, Wen Sun, Beifang Ning, Shanna Huang, Kun Wu, Chuanshu Huang, Mengchao Wu, Weifen Xie, and Hongyang Wang
Cancer Epidemiol. Biomarkers Prev., Jun 2009; 18: 1720 - 1729.

GAPDH depletion induces cell cycle arrest and resistance to antimetabolites in human carcinoma cell lines

Manali S Phadke, Natalia F Krynetskaia, Anurag K Mishra, and Evgeny Krynetskiy
J. Pharmacol. Exp. Ther., Jul 2009; 10.1124/jpet.109.155671.

A role for Chk1 in blocking transcriptional elongation of p21 RNA during the S-phase checkpoint

Rachel Beckerman, Aaron J. Donner, Melissa Mattia, Melissa J. Peart, James L. Manley, Joaquin M. Espinosa, and Carol Prives
Genes & Dev., Jun 2009; 23: 1364 - 1377.

HYPOXIA

Hypoxia-Inducible Factor 1 α Induces Fibrosis and Insulin Resistance in White Adipose Tissue

Nils Halberg, Tayeba Khan, Maria E. Trujillo, Ingrid Wernstedt-Asterholm, Alan D. Attie, Shariq Sherwani, Zhao V. Wang, Shira Landskroner-Eiger, Sean Dineen, Ulysses J. Magalang, Rolf A. Brekken, and Philipp E. Scherer
Mol. Cell. Biol., Aug 2009; 29: 4467 - 4483.

The molecular basis for impaired hypoxia-induced VEGF expression in diabetic tissues

Hariharan Thangarajah, Dachun Yao, Edward I. Chang, Yubin Shi, Leila Jazayeri, Ivan N. Vial, Robert D. Galiano, Xue-Liang Du, Raymon Grogan, Michael G. Galvez, Michael Januszyk, Michael Brownlee, and Geoffrey C. Gurtner
PNAS, Aug 2009; 106: 13505 - 13510.

Selective Inhibition of Hypoxia-Inducible Factor (HIF) Prolyl-Hydroxylase 1 Mediates Neuroprotection against Normoxic Oxidative Death via HIF- and CREB-Independent Pathways

Ambreena Siddiq, Leila R. Aminova, Carol M. Troy, Kyungsun Suh, Zachary Messer, Gregg L. Semenza, and Rajiv R. Ratan
J. Neurosci., Jul 2009; 29: 8828 - 8838.

Stretch-stimulated glucose uptake in skeletal muscle is mediated by reactive oxygen species and p38 MAP-kinase

Melissa A. Chambers, Jennifer S. Moylan, Jeffrey D. Smith, Laurie J. Goodyear, and Michael B. Reid
J. Physiol., Jul 2009; 587: 3363 - 3373.

Elevated hydrostatic pressure activates sodium/hydrogen exchanger-1 in rat optic nerve head astrocytes

Amritlal Mandal, Mohammad Shahidullah, Nicholas A. Delamere, and Marcos A. Terán
Am J Physiol Cell Physiol, Jul 2009; 297: C111 - C120

ION CHANNELS

Transient Receptor Potential Canonical 5 Channels Activate Ca²⁺/Calmodulin Kinase γ to Promote Axon Formation in Hippocampal Neurons

Monika A. Davare, Dale A. Fortin, Takeo Saneyoshi, Sean Nygaard, Stefanie Kaeck, Gary Banker, Thomas R. Soderling, and Gary A. Wayman
J. Neurosci., Aug 2009; 29: 9794 - 9808.

ASIC2 Subunits Target Acid-Sensing Ion Channels to the Synapse via an Association with PSD-95

Xiang-ming Zha, Vivian Costa, Anne Marie S. Harding, Leah Reznikov, Christopher J. Benson, and Michael J. Welsh
J. Neurosci., Jul 2009; 29: 8438 - 8446.

Phospholipase C- γ Binds Directly to the Na⁺/H⁺ Exchanger 3 and Is Required for Calcium Regulation of Exchange Activity

Nicholas C. Zachos, Damian B. van Rossum, Xuhang Li, Gabriela Caraveo, Rafiquel Sarker, Boyoung Cha, Sachin Mohan, Stephen Desiderio, Randen L. Patterson, and Mark Donowitz
J. Biol. Chem., Jul 2009; 284: 19437 - 19444.

METABOLISM

Widespread reorganization of metabolic enzymes into reversible assemblies upon nutrient starvation

Rammohan Narayanaswamy, Matthew Levy, Mark Tsechansky, Gwendolyn M. Stovall, Jeremy D. O'Connell, Jennifer Mirrielees, Andrew D. Ellington, and Edward M. Marcotte
PNAS, Jun 2009; 106: 10147 - 10152.

Attenuated suppression of the oxidative burst by cells dying in the presence of oxidized low density lipoprotein

Dmitry Namgaladze, Carla Jennewein, Stefan Preiss, Andreas von Knethen, and Bernhard Brüne
J. Lipid Res., June 2009; 10.1194/jlr.M800615-JLR200.

Depot-specific effects of the PPAR agonist rosiglitazone on adipose tissue glucose uptake and metabolism

William T. Festuccia, Pierre-Gilles Blanchard, Véronique Turcotte, Mathieu Laplante, Meltem Sariahmetoglu, David N. Brindley, and Yves Deshaies
J. Lipid Res., Jun 2009; 50: 1185 - 1194.

MATRIX METALOPROTEASES**Osteoclast-Derived Matrix Metalloproteinase-7, but Not Matrix Metalloproteinase-9, Contributes to Tumor-Induced Osteolysis**

Sophie Thiolloy, Jennifer Halpern, Ginger E. Holt, Herbert S. Schwartz, Gregory R. Mundy, Lynn M. Matrisian, and Conor C. Lynch
Cancer Res., Aug 2009; 69: 6747 - 6755.

MUSCLE**TRPC1 and STIM1 mediate capacitative Ca²⁺ entry in mouse pulmonary arterial smooth muscle cells**

Lih Chyuan Ng, Mary D. McCormack, Judith A. Airey, Cherie A. Singer, Phillip S. Keller, Xiao-Ming Shen, and Joseph R. Hume
J. Physiol., Jun 2009; 587: 2429 - 2442.

Taurine supplementation increases skeletal muscle force production and protects muscle function during and after high-frequency *in vitro* stimulation

Craig A. Goodman, Deanna Horvath, Christos Stathis, Trevor Mori, Kevin Croft, Robyn M. Murphy, and Alan Hayes
J Appl Physiol, Jul 2009; 107: 144 - 154.

Stretch-stimulated glucose uptake in skeletal muscle is mediated by reactive oxygen species and p38 MAP-kinase

Melissa A. Chambers, Jennifer S. Moylan, Jeffrey D. Smith, Laurie J. Goodyear, and Michael B. Reid
J. Physiol., Jul 2009; 587: 3363 - 3373.

 γ -Actin is required for cytoskeletal maintenance but not development

Inna A. Belyantseva, Benjamin J. Perrin, Kevin J. Sonnemann, Mei Zhu, Ruben Stepanyan, JoAnn McGee, Gregory I. Frolenkov, Edward J. Walsh, Karen H. Friderici, Thomas B. Friedman, and James M. Ervasti
PNAS, Jun 2009; 106: 9703 - 9708.

NEUROSCIENCE**Nucleocytoplasmic Shuttling of the Adapter Protein SH2B1 β (SH2-B β) Is Required for Nerve Growth Factor (NGF)-Dependent Neurite Outgrowth and Enhancement of Expression of a Subset of NGF-Responsive Genes**

Travis J. Maures, Linyi Chen, and Christin Carter-Su
Mol. Endocrinol., Jul 2009; 23: 1077 - 1091.

5-HT₄ Receptor-Mediated Neuroprotection and Neurogenesis in the Enteric Nervous System of Adult Mice

Min-Tsai Liu, Yung-Hui Kuan, Jingwen Wang, René Hen, and Michael D. Gershon
J. Neurosci., Aug 2009; 29: 9683 - 9699.

Tumor Necrosis Factor-Receptor-associated Factor-4 Is a Positive Regulator of Transforming Growth Factor- β Signaling That Affects Neural Crest Formation

Tuzer Kalkan, Yasuno Iwasaki, Chong Yon Park, and Gerald H. Thomsen
Mol. Biol. Cell, Jul 2009; 20: 3436 - 3450.

Decreased Brain-Derived Neurotrophic Factor Depends on Amyloid Aggregation State in Transgenic Mouse Models of Alzheimer's Disease

Shiyong Peng, Diego J. Garzon, Monica Marchese, William Klein, Stephen D. Ginsberg, Beverly M. Francis, Howard T. J. Mount, Elliott J. Mufson, Ahmad Salehi, and Margaret Fahnstock
J. Neurosci., Jul 2009; 29: 9321 - 9329.

Biological Activity of Nerve Growth Factor Precursor Is Dependent upon Relative Levels of Its Receptors

Raheleh Masoudi, Maria S. Ioannou, Michael D. Coughlin, Promila Pagadala, Kenneth E. Neet, Oliver Clewes, Shelley J. Allen, David Dawbarn, and Margaret Fahnestock
J. Biol. Chem., Jul 2009; 284: 18424 - 18433.

Conditional Forebrain Inactivation of Nicastrin Causes Progressive Memory Impairment and Age-Related Neurodegeneration

Katsuhiko Tabuchi, Guiquan Chen, Thomas C. Südhof, and Jie Shen
J. Neurosci., Jun 2009; 29: 7290 - 7301.

CK2 negatively regulates $G\alpha_s$ signaling

Heike Rebholz, Akinori Nishi, Sabine Liebscher, Angus C. Nairn, Marc Flajolet, and Paul Greengard
PNAS, Aug 2009; 106: 14096 - 14101.

TNF- α Preconditioning Protects Neurons via Neuron-Specific Up-Regulation of CREB-Binding Protein

Ramendra N. Saha, Anamitra Ghosh, Carlos A. Palencia, Yiu K. Fung, Serena M. Dudek, and Kalipada Pahan
J. Immunol., Aug 2009; 183: 2068 - 2078.

Mediobasal Hypothalamic Leucine Sensing Regulates Food Intake through Activation of a Hypothalamus–Brainstem Circuit

Clémence Blouet, Young-Hwan Jo, Xiaosong Li, and Gary J. Schwartz
J. Neurosci., Jul 2009; 29: 8302 - 8311.

Tissue Profiling of the Mammalian Central Nervous System Using Human Antibody-based Proteomics

Jan Mulder, Erik Björling, Kalle Jonasson, Henrik Wernérus, Sophia Hober, Tomas Hökfelt, and Mathias Uhlén
Mol. Cell. Proteomics, Jul 2009; 8: 1612 - 1622.

NEPHROLOGY

Local modulation of the natriuretic peptide system in the rat remnant kidney

Carla Santos-Araújo, Roberto Roncon-Albuquerque, Jr, Mónica Moreira-Rodrigues, Tiago Henriques-Coelho, Janete Quelhas-Santos, Bernardo Faria, Benedita Sampaio-Maia, Adelino F. Leite-Moreira, and Manuel Pestana
Nephrol. Dial. Transplant., Jun 2009; 24: 1774 - 1782.

Connexin 30 Deficiency Impairs Renal Tubular ATP Release and Pressure Natriuresis

Arnold Sipos, Sarah L. Vargas, Ildikó Toma, Fiona Hanner, Klaus Willecke, and János Peti-Peterdi
J. Am. Soc. Nephrol., Aug 2009; 20: 1724 - 1732.

PHOSPHATASES

Activation of Src by Protein Tyrosine Phosphatase 1B Is Required for ErbB2 Transformation of Human Breast Epithelial Cells

Luis E. Arias-Romero, Sayanti Saha, Olga Villamar-Cruz, Shu-Chin Yip, Stephen P. Ethier, Zhong-Yin Zhang, and Jonathan Chernoff
Cancer Res., Jun 2009; 69: 4582 - 4588.

RECEPTORS

Long-range activation of FKBP51 transcription by the androgen receptor via distal intronic enhancers

Harri Makkonen, Miia Kauhanen, Ville Paakinaho, Tiina Jääskeläinen, and Jorma J. Palvimö
Nucleic Acids Res., Jul 2009; 37: 4135 - 4148.

Role of Molecular Chaperones in G Protein $\beta 5$ /Regulator of G Protein Signaling Dimer Assembly and G Protein β Dimer Specificity

Alyson C. Howlett, Amy J. Gray, Jesse M. Hunter, and Barry M. Willardson
J. Biol. Chem., Jun 2009; 284: 16386 - 16399.

Endosomal Endothelin-converting Enzyme-1: A REGULATOR OF β -ARRESTIN-DEPENDENT ERK SIGNALING

Graeme S. Cottrell, Benjamin E. Padilla, Silvia Amadesi, Daniel P. Poole, Jane E. Murphy, Markus Hardt, Dirk Roosterman, Martin Steinhoff, and Nigel W. Bunnett
J. Biol. Chem., AUG 2009; 284: 22411 - 22425.

Regulation of Platelet-derived Growth Factor Receptor Function by Integrin-associated Cell Surface Transglutaminase

Evgeny A. Zemskov, Elena Loukinova, Irina Mikhailenko, Richard A. Coleman, Dudley K. Strickland, and Alexey M. Belkin
J. Biol. Chem., Jun 2009; 284: 16693 - 16703.

RNAi**CBP-mediated acetylation of histone H3 lysine 27 antagonizes Drosophila Polycomb silencing**

Feng Tie, Rakhee Banerjee, Carl A. Stratton, Jayashree Prasad-Sinha, Vincent Stepanik, Andrei Zlobin, Manuel O. Diaz, Peter C. Scacheri, and Peter J. Harte
Development, Sep 2009; 136: 3131 - 3141.

Up-regulation of miR-21 by HER2/neu Signaling Promotes Cell Invasion

Tzu-Hsuan Huang, Fangting Wu, Gabriel B. Loeb, Ruby Hsu, Amy Heidersbach, Allison Brincat, Dai Horiuchi, Robert J. Lebbink, Yin-Yuan Mo, Andrei Goga, and Michael T. McManus
J. Biol. Chem., Jul 2009; 284: 18515 - 18524.

The human Dcn1-like protein DCNL3 promotes Cul3 neddylation at membranes

Nathalie Meyer-Schaller, Yang-Chieh Chou, Izabela Sumara, Dale D. O. Martin, Thimo Kurz, Nadja Katheder, Kay Hofmann, Luc G. Berthiaume, Frank Sicheri, and Matthias Peter
PNAS, Jul 2009; 106: 12365 - 12370.

MicroRNA-125a-5p partly regulates the inflammatory response, lipid uptake, and ORP9 expression in oxLDL-stimulated monocyte/macrophages

Ting Chen, Zhouqing Huang, Liansheng Wang, Yue Wang, Feizhen Wu, Shu Meng, and Changqian Wang
Cardiovasc Res, Jul 2009; 83: 131 - 139.

MicroRNA Expression in Human Airway Smooth Muscle Cells: Role of miR-25 in Regulation of Airway Smooth Muscle Phenotype

Andrew R Kuhn, Karen Schlauch, Ronna Lao, Andrew J Halayko, William T Gerthoffer, and Cherie A Singer
Am. J. Respir. Cell Mol. Biol., Jun 2009; 10.1165/rcmb.2009-0123OC.

STEM CELLS**DNER, an Epigenetically Modulated Gene, Regulates Glioblastoma-Derived Neurosphere Cell Differentiation and Tumor Propagation**

Peng Sun, Shuli Xia, Bachchu Lal, Charles G. Eberhart, Alfredo Quinones-Hinojosa, Jarek Maciaczyk, William Matsui, Francesco DiMeco, Sara M. Piccirillo, Angelo L. Vescovi, and John Laterra
Stem Cells, Jul 2009; 27: 1473 - 1486.

Thrombin-cleaved osteopontin regulates hemopoietic stem and progenitor cell functions through interactions with $\alpha_9\beta_1$ and $\alpha_4\beta_1$ integrins

Jochen Grassinger, David N. Haylock, Melonie J. Storan, Gemma O. Haines, Brenda Williams, Genevieve A. Whitty, Andrew R. Vinson, Cheang Ly Be, Songhui Li, Esben S. Sørensen, Patrick P.L. Tam, David T. Denhardt, Dean Sheppard, Peter F. Choong, and Susan K. Nilsson
Blood, Jul 2009; 114: 49 - 59.

FOXO Transcription Factors Enforce Cell Cycle Checkpoints and Promote Survival of Hematopoietic Cells after DNA Damage

Hong Lei and Frederick W. Quelle
Mol. Cancer Res., Aug 2009; 7: 1294 - 1303

TRANSCRIPTION FACTORS

Long-range activation of FKBP51 transcription by the androgen receptor via distal intronic enhancers

Harri Makkonen, Miia Kauhanen, Ville Paakinaho, Tiina Jääskeläinen, and Jorma J. Palvimo
Nucleic Acids Res., Jul 2009; 37: 4135 - 4148.

TNF- α Preconditioning Protects Neurons via Neuron-Specific Up-Regulation of CREB-Binding Protein

Ramendra N. Saha, Anamitra Ghosh, Carlos A. Palencia, Yiu K. Fung, Serena M. Dudek, and Kalipada Pahan
J. Immunol., Aug 2009; 183: 2068 - 2078.

Role of GRP78/BiP Degradation and ER Stress in Deoxynivalenol-Induced Interleukin-6 Upregulation in the Macrophage

Yuhui Shi, Katie Porter, Narayanan Parameswaran, Hee Kyong Bae, and James J. Pestka
Toxicol. Sci., Jun 2009; 109: 247 - 255.

FOXO Transcription Factors Enforce Cell Cycle Checkpoints and Promote Survival of Hematopoietic Cells after DNA Damage

Hong Lei and Frederick W. Quelle
Mol. Cancer Res., Aug 2009; 7: 1294 – 1303

Insulin acts through FOXO3a to activate transcription of Plasminogen Activator Inhibitor Type 1

Ushma R. Jag, Jiri Zavadil, and Frederick M. Stanley
Mol. Endocrinol., Jul 2009; 10.1210/me.2008-0421.

Annexin-1 Regulates Macrophage IL-6 and TNF via Glucocorticoid-Induced Leucine Zipper

Yuan H. Yang, Daniel Aeberli, April Dacumos, Jin R. Xue, and Eric F. Morand
J. Immunol., Jul 2009; 183: 1435 - 1445.

UBIQUITINATION AND SUMOYLATION

Syk Tyrosine 317 Negatively Regulates Osteoclast Function via the Ubiquitin-Protein Isopeptide Ligase Activity of Cbl

Wei Zou, Jennifer L. Reeve, Haibo Zhao, F. Patrick Ross, and Steven L. Teitelbaum
J. Biol. Chem., Jul 2009; 284: 18833 - 18839.

Role of SUMO in RNF4-mediated Promyelocytic Leukemia Protein (PML) Degradation: Sumoylation of pml and phospho-switch control of its sumo binding domain dissected in living cells

Yann Percherancier, Delphine Germain-Desprez, Frédéric Galisson, Xavier H. Mascle, Laurent Dianoux, Patricia Estephan, Mounira K. Chelbi-Alix, and Muriel Aubry
J. Biol. Chem., Jun 2009; 284: 16595 - 16608.

TRIM32 is an E3 ubiquitin ligase for dysbindin

Matthew Locke, Caroline L. Tinsley, Matthew A. Benson, and Derek J. Blake
Hum. Mol. Genet., Jul 2009; 18: 2344 - 2358.

Degradation of regulator of calcineurin 1 (RCAN1) is mediated by both chaperone-mediated autophagy and ubiquitin proteasome pathways

Heng Liu, Pin Wang, Weihong Song, and Xiulian Sun
FASEB J, Jun 2009; 10.1096/fj.09-134296.

The human Dcn1-like protein DCNL3 promotes Cul3 neddylation at membranes

Nathalie Meyer-Schaller, Yang-Chieh Chou, Izabela Sumara, Dale D. O. Martin, Thimo Kurz, Nadja Katheder, Kay Hofmann, Luc G. Berthiaume, Frank Sicheri, and Matthias Peter
PNAS, Jul 2009; 106: 12365 - 12370.

The fusion proteins TEL-PDGFR β and FIP1L1-PDGFR α escape ubiquitination and degradation

Federica Toffalini, Anders Kallin, Peter Vandenberghe, Pascal Pierre, Lucienne Michaux, Jan Cools, and Jean-Baptiste Demoulin
Haematologica, Aug 2009; 94: 1085 - 1093.

VIROLOGY

Murine Cytomegalovirus Capsid Assembly Is Dependent on US22 Family Gene M140 in Infected Macrophages

Laura K. Hanson, Jacquelyn S. Slater, Victoria J. Cavanaugh, William W. Newcomb, Lisa L. Bolin, Christine N. Nelson, Lisa D. Fetters, Qiyi Tang, Jay C. Brown, Gerd G. Maul, and Ann E. Campbell
J. Virol., Aug 2009; 83: 7449 - 7456.

Clathrin adaptor AP1B controls adenovirus infectivity of epithelial cells

Fernando Diaz, Diego Gravotta, Ami Deora, Ryan Schreiner, John Schoggins, Erik Falck-Pedersen, and Enrique Rodriguez-Boulan
PNAS, Jul 2009; 106: 11143 - 11148.

Vaccinia Virus H7 Protein Contributes to the Formation of Crescent Membrane Precursors of Immature Virions

P. S. Satheshkumar, Andrea Weisberg, and Bernard Moss
J. Virol., Sep 2009; 83: 8439 - 8450.

IL-27 inhibition of HIV-1 involves an intermediate induction of type I IFN

Teresa Greenwell-Wild, Nancy Vazquez, Wenwen Jin, Zoila Rangel, Peter Munson, and Sharon M. Wahl
Blood, Jun 2009; 10.1182/blood-2009-03-211540.

Matrix Mediates the Functional Link between Human Immunodeficiency Virus Type 1 RNA Nuclear Export Elements and the Assembly Competency of Gag in Murine Cells

Nathan M. Sherer, Chad M. Swanson, Stelios Papaioannou, and Michael H. Malim
J. Virol., Sep 2009; 83: 8525 - 8535

Ebola Virus VP35 Antagonizes PKR Activity through Its C-Terminal Interferon Inhibitory Domain

Michael Schümann, Thorsten Gantke, and Elke Mühlberger
J. Virol., Sep 2009; 83: 8993 - 8997.

Latency-Associated Nuclear Antigen of Kaposi's Sarcoma-Associated Herpesvirus (KSHV) Upregulates Survivin Expression in KSHV-Associated B-Lymphoma Cells and Contributes to Their Proliferation

Jie Lu, Subhash C. Verma, Masanao Murakami, Qiliang Cai, Pankaj Kumar, Bingyi Xiao, and Erle S. Robertson
J. Virol., Jul 2009; 83: 7129 - 7141.

Human Papillomavirus Type 16 Infection of Human Keratinocytes Requires Clathrin and Caveolin-1 and Is Brefeldin A Sensitive

Valerie Laniosz, Sarah A. Dabydeen, Mallory A. Havens, and Patricio I. Meneses
J. Virol., Aug 2009; 83: 8221 - 8232.

Potent neutralization of anthrax edema toxin by a humanized monoclonal antibody that competes with calmodulin for edema factor binding

Zhaochun Chen, Mahtab Moayeri, Huaying Zhao, Devorah Crown, Stephen H. Leppla, and Robert H. Purcell
PNAS, Aug 2009; 106: 13487 - 13492.

Proanthocyanidin from Blueberry Leaves Suppresses Expression of Subgenomic Hepatitis C Virus RNA

Masahiko Takeshita, Yo-ichi Ishida, Ena Akamatsu, Yusuke Ohmori, Masayuki Sudoh, Hirofumi Uto, Hirohito Tsubouchi, and Hiroaki Kataoka
J. Biol. Chem., Aug 2009; 284: 21165 - 21176.

APPLICATIONS

In-Cell Western

Activation of Src by Protein Tyrosine Phosphatase 1B Is Required for ErbB2 Transformation of Human Breast Epithelial Cells

Luis E. Arias-Romero, Sayanti Saha, Olga Villamar-Cruz, Shu-Chin Yip, Stephen P. Ethier, Zhong-Yin Zhang, and Jonathan Chernoff
Cancer Res., Jun 2009; 69: 4582 - 4588.

Pathogenic role of angiotensin II and oxidized LDL in obstructive sleep apnoea

T. Kizawa, Y. Nakamura, S. Takahashi, S. Sakurai, K. Yamauchi, and H. Inoue
Eur. Respir. J., Jul 2009; 10.1183/09031936.00009709.

Coupling of Ionic Events to Protein Kinase Signaling Cascades upon Activation of 7 Nicotinic Receptor: COOPERATIVE REGULATION OF α_2 -INTEGRIN EXPRESSION AND Rho KINASE ACTIVITY

Alexander I. Chernyavsky, Juan Arredondo, Jing Qian, Valentin Galitovskiy, and Sergei A. Grando
J. Biol. Chem., Aug 2009; 284: 22140 - 22148.

Agonist-biased signaling via Proteinase Activated Receptor-2: differential activation of calcium and MAPkinase pathways.

Rithwik Ramachandran, Koichiro Mihara, Maneesh Mathur, Moulay Driss Rochdi, Michel Bouvier, Kathryn DeFea, and Morley D. Hollenberg
Mol. Pharmacol., Jul 2009; 10.1124/mol.109.055509.

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)

Quantitative Western (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

MousePod

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)

Quantitative Reverse Phase Protein Arrays

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.

DNA Methylation Assay

Quantification of global DNA Methylation with infrared fluorescence in liver and muscle tissues of differentially fed boars

H. Braunschweig
Luminescence 2009 (www.interscience.wiley.com) DOI 10.1002/bio.1098)

LI-COR and Odyssey are registered trademarks of LI-COR, Inc.
All other trademarks belong to their respective owners.



4647 Superior Street • P.O. Box 4000 • Lincoln, Nebraska 68504 USA
North America: 800-645-4267 • International: 402-467-0700 • FAX: 402-467-0819
LI-COR GmbH Germany, Serving Europe and Africa: +49 (0) 6172 17 17 771
LI-COR UK Ltd. UK, Serving UK, Ireland, Scandinavia: +44 (0) 1223 422104
www.licor.com

Doc #979-07647
Rev 0909