

1: Diggin' on U(biquitin): A Novel Method for the Identification of Physiological E3 Ubiquitin Ligase Substrates

Carrie E. Rubel, Jonathan C. Schisler, Eric D. Hamlett, Robert M. DeKroon, Mathias Gautel, Oscar Alzate, Cam Patterson

Cell Biochem Biophys. Author manuscript; available in PMC 2014 September 1. Published in final edited form as: Cell Biochem Biophys. 2013 September; 67(1): 127–138.

doi: 10.1007/s12013-013-9624-6

PMCID: PMC3758785

2: Recombinant Scorpine Produced Using SUMO Fusion Partner in Escherichia coli Has the Activities against Clinically Isolated Bacteria and Inhibits the Plasmodium falciparum Parasitemia In Vitro

Chao Zhang, Xinlong He, Yaping Gu, Huayun Zhou, Jun Cao, Qi Gao

PLoS One. 2014; 9(7): e103456. Published online 2014 July 28.

doi: 10.1371/journal.pone.0103456

PMCID: PMC4113386

3: The Aspartate-Less Receiver (ALR) Domains: Distribution, Structure and Function

Andrew F. Maule, David P. Wright, Joshua J. Weiner, Lanlan Han, Francis C. Peterson, Brian F. Volkman, Nicholas R. Silvaggi, Andrew T. Ulijasz

PLoS Pathog. 2015 April; 11(4): e1004795. Published online 2015 April 13.

doi: 10.1371/journal.ppat.1004795 Correction in: PLoS Pathog. 2015 June; 11(6): e1004951.

PMCID: PMC4395418

4: BLAP-Tags, TUBEs and DUB-Chips: Combined Novel Technologies will Advance Molecular Epithelial Physiology

Kirk L. Hamilton

Front Physiol. 2012; 3: 137. Published online 2012 May 15. doi: 10.3389/fphys.2012.00137

PMCID: PMC3351676

5: A Comparative Analysis of the Ubiquitination Kinetics of Multiple Degrons to Identify an Ideal Targeting Sequence for a Proteasome Reporter

Adam T. Melvin, Gregory S. Woss, Jessica H. Park, Lukas D. Dumberger, Marcey L. Waters, Nancy L. Allbritton

PLoS One. 2013; 8(10): e78082. Published online 2013 October 29.

doi: 10.1371/journal.pone.0078082

PMCID: PMC3812159

6: The LRR and RING Domain Protein LRSAM1 Is an E3 Ligase Crucial for Ubiquitin-Dependent Autophagy of Intracellular Salmonella Typhimurium

Alan Huett, Robert J. Heath, Jakob Begun, Slim O. Sassi, Leigh A. Baxt, Jatin M. Vyas, Marcia B. Goldberg, Ramnik J. Xavier

Cell Host Microbe. Author manuscript; available in PMC 2013 December 13. Published in final edited form as: Cell Host Microbe. 2012 December 13; 12(6): 778–790.

doi: 10.1016/j.chom.2012.10.019

PMCID: PMC3785244

- 7: Methods for Quantification of in vivo Changes in Protein Ubiquitination following Proteasome and Deubiquitinase Inhibition
Namrata D. Udeshi, D. R. Mani, Thomas Eisenhaure, Philipp Mertins, Jacob D. Jaffe, Karl R. Clauser, Nir Hacohen, Steven A. Carr
Mol Cell Proteomics. 2012 May; 11(5): 148–159. Published online 2012 April 14.
doi: 10.1074/mcp.M111.016857
PMCID: PMC3418844
- 8: MDM2 Mediates Nonproteolytic Polyubiquitylation of the DEAD-Box RNA Helicase DDX24
Takayoshi Yamauchi, Masaaki Nishiyama, Toshiro Moroishi, Kanae Yumimoto, Keiichi I. Nakayama
Mol Cell Biol. 2014 September; 34(17): 3321–3340. doi: 10.1128/MCB.00320-14
PMCID: PMC4135551
- 9: Deubiquitylase, DeSUMOylase, and DeISGylase Activity Microarrays for Assay of Substrate Preference and Functional Modifiers
Christian M. Loch, Charles L. Cuccherini, Craig A. Leach, James E. Strickler
Mol Cell Proteomics. 2011 January; 10(1): M110.002402. Published online 2010 October 18.
doi: 10.1074/mcp.M110.002402
PMCID: PMC3013450
- 10: Facilitated Tau Degradation by USP14 Aptamers via Enhanced Proteasome Activity
Jung Hoon Lee, Seung Kyun Shin, Yanxialei Jiang, Won Hoon Choi, Chaesun Hong, Dong-Eun Kim, Min Jae Lee
Sci Rep. 2015; 5: 10757. Published online 2015 June 4. doi: 10.1038/srep10757
PMCID: PMC4455164
- 11: Role of ubiquitylation and USP8-dependent deubiquitylation in the endocytosis and lysosomal targeting of plasma membrane KCa3.1
Corina M. Balut, Christian M. Loch, Daniel C. Devor
FASEB J. 2011 November; 25(11): 3938–3948. doi: 10.1096/fj.11-187005
PMCID: PMC3205838
- 12: Omega 3 fatty acids chemosensitize multidrug resistant colon cancer cells by down-regulating cholesterol synthesis and altering detergent resistant membranes composition
Giada Gelsomino, Paola A Corsetto, Ivana Campia, Gigliola Montorfano, Joanna Kopecka, Barbara Castella, Elena Gazzano, Dario Ghigo, Angela M Rizzo, Chiara Riganti
Mol Cancer. 2013; 12: 137. Published online 2013 November 13. doi: 10.1186/1476-4598-12-137
PMCID: PMC4225767
- 13: Dissecting the interactions of SERRATE with RNA and DICER-LIKE 1 in Arabidopsis microRNA precursor processing
Yuji Iwata, Masateru Takahashi, Nina V. Fedoroff, Samir M. Hamdan

Nucleic Acids Res. 2013 October; 41(19): 9129–9140. Published online 2013 August 5.

doi: 10.1093/nar/gkt667

PMCID: PMC3799435

14: Factor Xa Activation of Factor V is of Paramount Importance in Initiating the Coagulation System: Lessons from a Tick Salivary Protein

Tim J. Schuijt, Kamran Bakhtiari, Sirlei Daffre, Kathleen DePonte, Simone J.H. Wielders, J. Arnoud Marquart, Joppe W. Hovius, Tom van der Poll, Erol Fikrig, Matthew W. Bunce, Rodney M. Camire, Gerry A.F. Nicolaes, Joost C.M. Meijers, Cornelis van 't Veer

Circulation. Author manuscript; available in PMC 2014 July 16. Published in final edited form as: Circulation. 2013 July 16; 128(3): 10.1161/CIRCULATIONAHA.113.003191. Published online 2013 July 1. doi: 10.1161/CIRCULATIONAHA.113.003191

PMCID: PMC3826089

15: Structural and Functional Characterization of MppR, an Enduracididine Biosynthetic Enzyme from Streptomyces hygroscopicus: Functional Diversity in the Acetoacetate Decarboxylase-Like Superfamily

A. Maxwell Burroughs, Robert W. Hoppe, Neal C. Goebel, Bilal H. Sayyed, Tyler J. Voegtline, Alan W. Schwabacher, T. Mark Zabriskie, Nicholas R. Silvaggi

Biochemistry. Author manuscript; available in PMC 2014 July 2. Published in final edited form as: Biochemistry. 2013 July 2; 52(26): 4492–4506. Published online 2013 June 21.

doi: 10.1021/bi400397k

PMCID: PMC3743547

16: Chronic carvedilol treatment partially reverses the right ventricular failure transcriptional profile in experimental pulmonary hypertension

Jennifer I. Drake, Jose Gomez-Arroyo, Catherine I. Dumur, Donatas Kraskauskas, Ramesh Natarajan, Harm J. Bogaard, Paul Fawcett, Norbert F. Voelkel

Physiol Genomics. 2013 June 15; 45(12): 449–461. Published online 2013 April 30.

doi: 10.1152/physiolgenomics.00166.2012

PMCID: PMC3680777

17: A Co-Translational Ubiquitination Pathway For Quality Control of Misfolded Proteins
Feng Wang, Larissa A. Durfee, Jon M. Huibregtse

Mol Cell. Author manuscript; available in PMC 2014 May 9. Published in final edited form as: Mol Cell. 2013 May 9; 50(3): 368–378. Published online 2013 April 11.

doi: 10.1016/j.molcel.2013.03.009

PMCID: PMC3654026

18: MreA Functions in the Global Regulation of Methanogenic Pathways in Methanosarcina acetivorans

21: Identification of Targeting Peptides for Ischemic Myocardium by In Vivo Phage Display
Sachiko Kanki, Diana E. Jaalouk, Samuel Lee, Alvin Y.C. Yu, Joseph Gannon, Richard T. Lee

J Mol Cell Cardiol. Author manuscript; available in PMC 2012 May 1. Published in final edited form as: J Mol Cell Cardiol. 2011 May; 50(5): 841–848. Published online 2011 February 24.
doi: 10.1016/j.yjmcc.2011.02.003
PMCID: PMC3075368

22: Inhibition of UCH-L1 in oligodendroglial cells results in microtubule stabilization and prevents α -synuclein aggregate formation by activating the autophagic pathway: implications for multiple system atrophy
Katharina Pukaß, Christiane Richter-Landsberg
Front Cell Neurosci. 2015; 9: 163. Published online 2015 May 5. doi: 10.3389/fncel.2015.00163
PMCID: PMC4419839

23: Identification of a Tsetse Fly Salivary Protein with Dual Inhibitory Action on Human Platelet Aggregation
Guy Caljon, Karin De Ridder, Patrick De Baetselier, Marc Coosemans, Jan Van Den Abbeele
PLoS One. 2010; 5(3): e9671. Published online 2010 March 23.
doi: 10.1371/journal.pone.0009671
PMCID: PMC2843633

24: Biosimilars: the process is the product. The example of recombinant streptokinase
C Thelwell, C Longstaff
J Thromb Haemost. 2014 August; 12(8): 1229–1233. Published online 2014 July 31.
doi: 10.1111/jth.12629
PMCID: PMC4241032

25: Recombinant Antigens Expressed in *Pichia pastoris* for the Diagnosis of Sleeping Sickness Caused by *Trypanosoma brucei gambiense*
Stijn Rogé, Liesbeth Van Nieuwenhove, Magali Meul, Annick Heykers, Annette Brouwer de Koning, Nicolas Bebronne, Yves Guisez, Philippe Büscher
PLoS Negl Trop Dis. 2014 July; 8(7): e3006. Published online 2014 July 17.
doi: 10.1371/journal.pntd.0003006
PMCID: PMC4102443

26: Anterograde Trafficking of KCa3.1 in Polarized Epithelia Is Rab1- and Rab8-Dependent and Recycling Endosome-Independent
Claudia A. Bertuccio, Shih-Liang Lee, Guangyu Wu, Michael B. Butterworth, Kirk L. Hamilton, Daniel C. Devor
PLoS One. 2014; 9(3): e92013. Published online 2014 March 14.
doi: 10.1371/journal.pone.0092013
PMCID: PMC3954861

27: MDM2 restrains estrogen-mediated AKT activation by promoting TBK1-dependent HPIP degradation
K Shostak, F Patrascu, S I Göktuna, P Close, L Borgs, L Nguyen, F Olivier, A Rammal, H Brinkhaus, M Bentires-Alj, J-C Marine, A Chariot
Cell Death Differ. 2014 May; 21(5): 811–824. Published online 2014 January 31.
doi: 10.1038/cdd.2014.2

PMCID: PMC3978309

28: Ataxia and hypogonadism caused by the loss of ubiquitin ligase activity of the U box protein CHIP

Chang-He Shi, Jonathan C. Schisler, Carrie E. Rubel, Song Tan, Bo Song, Holly McDonough, Lei Xu, Andrea L. Portbury, Cheng-Yuan Mao, Cadence True, Rui-Hao Wang, Qing-Zhi Wang, Shi-Lei Sun, Stephanie B. Seminara, Cam Patterson, Yu-Ming Xu

Hum Mol Genet. 2014 February 15; 23(4): 1013–1024. Published online 2013 October 9.

doi: 10.1093/hmg/ddt497

PMCID: PMC3900109

29: Restoring the Procofactor State of Factor Va-like Variants by Complementation with B-domain Peptides

Matthew W. Bunce, Mettine H. A. Bos, Sriram Krishnaswamy, Rodney M. Camire

J Biol Chem. 2013 October 18; 288(42): 30151–30160. Published online 2013 September 6.

doi: 10.1074/jbc.M113.506840

PMCID: PMC3798483

30: To fuse or not to fuse: What is your purpose?

Mark R Bell, Mark J Engleka, Asim Malik, James E Strickler

Protein Sci. 2013 November; 22(11): 1466–1477. Published online 2013 August 26.

doi: 10.1002/pro.2356

PMCID: PMC3831663

31: S-nitrosylation regulates mitochondrial quality control via activation of parkin

Kentaro Ozawa, Akira T. Komatsubara, Yuhei Nishimura, Tomoyo Sawada, Hiroto Kawafune, Hiroki Tsumoto, Yuichi Tsuji, Jing Zhao, Yoji Kyotani, Toshio Tanaka, Ryosuke Takahashi, Masanori Yoshizumi

Sci Rep. 2013; 3: 2202. Published online 2013 July 16. doi: 10.1038/srep02202

PMCID: PMC3712319

32: Purification, crystallization and preliminary X-ray analysis of TP0435 (Tp17) from the syphilis spirochete *Treponema pallidum*

Chad A. Brautigam, Ranjit K. Deka, Michael V. Norgard

Acta Crystallogr Sect F Struct Biol Cryst Commun. 2013 April 1; 69(Pt 4): 453–455. Published online 2013 March 28. doi: 10.1107/S1744309113006246

PMCID: PMC3614177

33: The TP0796 Lipoprotein of *Treponema pallidum* Is a Bimetal-dependent FAD Pyrophosphatase with a Potential Role in Flavin Homeostasis

Ranjit K. Deka, Chad A. Brautigam, Wei Z. Liu, Diana R. Tomchick, Michael V. Norgard

J Biol Chem. 2013 April 19; 288(16): 11106–11121. Published online 2013 February 27.

doi: 10.1074/jbc.M113.449975

PMCID: PMC3630870

34: Solution Structure of Escherichia coli FeoA and Its Potential Role in Bacterial Ferrous Iron Transport

Cheryl K. Y. Lau, Hiroaki Ishida, Zhihong Liu, Hans J. Vogel
J Bacteriol. 2013 January; 195(1): 46–55. doi: 10.1128/JB.01121-12
PMCID: PMC3536175

35: Enhanced protein expression in mammalian cells using engineered SUMO fusions: Secreted phospholipase A2

Raymond J. Peroutka, Nabil Elshourbagy, Tara Piech, Tauseef R. Butt
Protein Sci. 2008 September; 17(9): 1586–1595. doi: 10.1110/ps.035576.108
PMCID: PMC2525526

36: Bioluminescence Assay Platform for Selective and Sensitive Detection of Ub/Ubl Proteases

Steven J. Orcutt, Jian Wu, Michael J. Eddins, Craig A. Leach, James E. Strickler
Biochim Biophys Acta. Author manuscript; available in PMC 2013 November 1. Published in final edited form as: Biochim Biophys Acta. 2012 November; 1823(11): 2079–2086. Published online 2012 June 15. doi: 10.1016/j.bbamcr.2012.06.004
PMCID: PMC3465522

37: Structural, bioinformatic, and in vivo analyses of two Treponema pallidum lipoproteins reveal a unique TRAP transporter

Ranjit K. Deka, Chad A. Brautigam, Martin Goldberg, Peter Schuck, Diana R. Tomchick, Michael V. Norgard
J Mol Biol. Author manuscript; available in PMC 2013 March 9. Published in final edited form as: J Mol Biol. 2012 March 9; 416(5): 678–696. Published online 2012 January 27. doi: 10.1016/j.jmb.2012.01.015
PMCID: PMC3289903

38: Characterization of Selective Ubiquitin and Ubiquitin-Like Protease Inhibitors Using a Fluorescence-Based Multiplex Assay Format

Xufan Tian, Nigora S. Isamiddinova, Raymond J. Peroutka, Seth J. Goldenberg, Michael R. Mattern, Benjamin Nicholson, Craig Leach
Assay Drug Dev Technol. 2011 April; 9(2): 165–173. doi: 10.1089/adt.2010.0317
PMCID: PMC3065724

39: Efficient and rapid protein expression and purification of small high disulfide containing sweet protein brazzein in E. coli

Fariba Assadi-Porter, Sammy Patry, John L. Markley
Protein Expr Purif. Author manuscript; available in PMC 2009 April 1. Published in final edited form as: Protein Expr Purif. 2008 April; 58(2): 263–268. Published online 2007 December 3. doi: 10.1016/j.pep.2007.11.009
PMCID: PMC2374762

40: Evidence that Altered Cis Element Spacing Affects PpsR Mediated Redox Control of Photosynthesis Gene Expression in Rubrivivax gelatinosus

Takayuki Shimizu, Zhuo Cheng, Katsumi Matsuura, Shinji Masuda, Carl E. Bauer

PLoS One. 2015; 10(6): e0128446. Published online 2015 June 1.
doi: 10.1371/journal.pone.0128446
PMCID: PMC4452267

41: Development of two novel high-throughput assays to quantify ubiquitylated proteins in cell lysates: application to screening of new anti-malarials
Lydia Mata-Cantero, Concepción Cid, Maria G Gomez-Lorenzo, Wendy Xolalpa, Fabienne Aillet, J Julio Martín, Manuel S Rodriguez
Malar J. 2015; 14: 200. Published online 2015 May 14. doi: 10.1186/s12936-015-0708-1
PMCID: PMC4440562

42: A secretory kinase complex regulates extracellular protein phosphorylation
Jixin Cui, Junyu Xiao, Vincent S Tagliabracci, Jianzhong Wen, Meghdad Rahdar, Jack E Dixon
eLife. 2015; 4: e06120. Published online 2015 March 19. doi: 10.7554/eLife.06120Includes additional comments & authors
PMCID: PMC4421793

43: Lesion bypass activity of DNA polymerase θ (POLQ) is an intrinsic property of the pol domain and depends on unique sequence inserts
Matthew Hogg, Mineaki Seki, Richard D. Wood, Sylvie Doublé, Susan S. Wallace
J Mol Biol. Author manuscript; available in PMC 2012 January 21. Published in final edited form as: J Mol Biol. 2011 January 21; 405(3): 642–652. Published online 2010 November 2.
doi: 10.1016/j.jmb.2010.10.041Correction in: volume 412 on page 316
PMCID: PMC3025778

44: HIV-1 Vpr Loads Uracil DNA Glycosylase-2 onto DCAF1, a Substrate Recognition Subunit of a Cullin 4A-RING E3 Ubiquitin Ligase for Proteasome-dependent Degradation
Jinwoo Ahn, Thomas Vu, Zach Novince, Jennifer Guerrero-Santoro, Vesna Rapic-Otrin, Angela M. Gronenborn
J Biol Chem. 2010 November 26; 285(48): 37333–37341. Published online 2010 September 24.
doi: 10.1074/jbc.M110.133181
PMCID: PMC2988339

45: A novel RNA molecular signature for activation of 2'-5' oligoadenylate synthetase-1
Virginia K. Vachon, Brenda M. Calderon, Graeme L. Conn
Nucleic Acids Res. 2015 January 9; 43(1): 544–552. Published online 2014 December 4.
doi: 10.1093/nar/gku1289
PMCID: PMC4288181

46: Ubiquitin-proteasome system controls ciliogenesis at the initial step of axoneme extension
Kousuke Kasahara, Yoshitaka Kawakami, Tohru Kiyono, Shigenobu Yonemura, Yoshifumi Kawamura, Saho Era, Fumio Matsuzaki, Naoki Goshima, Masaki Inagaki
Nat Commun. 2014 October 1; 5: 5081. Published online 2014 October 1.
doi: 10.1038/ncomms6081
PMCID: PMC4205846

47: Solution Structure of the 2A Protease from a Common Cold Agent, Human Rhinovirus C2, Strain W12

Woonghee Lee, Kelly E. Watters, Andrew T. Troupis, Nichole M. Reinen, Fabian P. Suchy, Kylie L. Moyer, Ronnie O. Frederick, Marco Tonelli, David J. Aceti, Ann C. Palmenberg, John L. Markley

PLoS One. 2014; 9(6): e97198. Published online 2014 June 17.

doi: 10.1371/journal.pone.0097198

PMCID: PMC4061012

48: Nonequilibrium dynamics and ultraslow relaxation of confined DNA during viral packaging

Zachary T. Berndsen, Nicholas Keller, Shelley Grimes, Paul J. Jardine, Douglas E. Smith

Proc Natl Acad Sci U S A. 2014 June 10; 111(23): 8345–8350. Published online 2014 May 27.

doi: 10.1073/pnas.1405109111

PMCID: PMC4060647

49: Membrane protein stability can be compromised by detergent interactions with the extramembranous soluble domains

Zhengrong Yang, Chi Wang, Qingxian Zhou, Jianli An, Ellen Hildebrandt, Luba A Aleksandrov, John C Kappes, Lawrence J DeLucas, John R Riordan, Ina L Urbatsch, John F Hunt, Christie G Brouillette

Protein Sci. 2014 June; 23(6): 769–789. Published online 2014 March 21. doi: 10.1002/pro.2460

PMCID: PMC4093953

50: Ubiquitin C-terminal Hydrolase L1 Interacts with Choline Transporter in Cholinergic Cells
Sigurd Hartnett, Fan Zhang, Allison Abitz, Yifan Li

Neurosci Lett. Author manuscript; available in PMC 2015 April 3. Published in final edited form as: Neurosci Lett. 2014 April 3; 564: 115–119. Published online 2014 February 11.

doi: 10.1016/j.neulet.2014.02.007

PMCID: PMC4024205

51: Structural and Computational Studies of the Staphylococcus aureus Sortase B-Substrate Complex Reveal a Substrate-stabilized Oxyanion Hole

Alex W. Jacobitz, Jeff Wereszczynski, Sung Wook Yi, Brendan R. Amer, Grace L. Huang, Angelyn V. Nguyen, Michael R. Sawaya, Michael E. Jung, J. Andrew McCammon, Robert T. Clubb

J Biol Chem. 2014 March 28; 289(13): 8891–8902. Published online 2014 February 11.

doi: 10.1074/jbc.M113.509273

PMCID: PMC3979406

52: Activation-triggered subunit exchange between CaMKII holoenzymes facilitates the spread of kinase activity

Margaret Stratton, Il-Hyung Lee, Moitrayee Bhattacharyya, Sune M Christensen, Luke H Chao, Howard Schulman, Jay T Groves, John Kuriyan

eLife. 2014; 3: e01610. Published online 2014 January 28. doi: 10.7554/eLife.01610 Includes additional comments & authors Correction in: eLife. 2014; 3: e02490.

PMCID: PMC3901001

53: Vitamin B12 regulates photosystem gene expression via the CrtJ antirepressor AerR in *Rhodobacter capsulatus*

Zhuo Cheng, Keran Li, Loubna A. Hammad, Jonathan A. Karty, Carl E. Bauer

Mol Microbiol. Author manuscript; available in PMC 2015 February 1. Published in final edited form as: Mol Microbiol. 2014 February; 91(4): 649–664. Published online 2014 January 9.

doi: 10.1111/mmi.12491

PMCID: PMC3946051

54: Differential Loss of Prolyl Isomerase or Chaperone Activity of Ran-binding Protein 2 (Ranbp2) Unveils Distinct Physiological Roles of Its Cyclophilin Domain in Proteostasis

Kyoung-in Cho, Hemangi Patil, Eugene Senda, Jessica Wang, Haiqing Yi, Sunny Qiu, Dosuk Yoon, Minzhong Yu, Andrew Orry, Neal S. Peachey, Paulo A. Ferreira

J Biol Chem. 2014 February 21; 289(8): 4600–4625. Published online 2014 January 8.

doi: 10.1074/jbc.M113.538215

PMCID: PMC3931022

55: Structure-guided Development of Specific Pyruvate Dehydrogenase Kinase Inhibitors Targeting the ATP-binding Pocket

Shih-Chia Tso, Xiangbing Qi, Wen-Jun Gui, Cheng-Yang Wu, Jacinta L. Chuang, Ingrid Wernstedt-Asterholm, Lorraine K. Morlock, Kyle R. Owens, Philipp E. Scherer, Noelle S. Williams, Uttam K. Tambar, R. Max Wynn, David T. Chuang

J Biol Chem. 2014 February 14; 289(7): 4432–4443. Published online 2013 December 19.

doi: 10.1074/jbc.M113.533885

PMCID: PMC3924305

56: Haploid Genetic Screens Identify an Essential Role for PLP2 in the Downregulation of Novel Plasma Membrane Targets by Viral E3 Ubiquitin Ligases

Richard T. Timms, Lidia M. Duncan, Iva A. Tchasonnikarova, Robin Antrobus, Duncan L. Smith, Gordon Dougan, Michael P. Weekes, Paul J. Lehner

PLoS Pathog. 2013 November; 9(11): e1003772. Published online 2013 November 21.

doi: 10.1371/journal.ppat.1003772

PMCID: PMC3836740

57: Substrate Binding Promotes Formation of the Skp1-Cul1-Fbx13 (SCFFbx13) Protein Complex

Kanae Yumimoto, Tetsuya Muneoka, Tomohiro Tsuboi, Keiichi I. Nakayama

J Biol Chem. 2013 November 8; 288(45): 32766–32776. Published online 2013 September 30.

doi: 10.1074/jbc.M113.511303

PMCID: PMC3820910

58: F-box and WD Repeat Domain-containing-7 (Fbxw7) Protein Targets Endoplasmic Reticulum-anchored Osteogenic and Chondrogenic Transcriptional Factors for Degradation

Kanae Yumimoto, Masaki Matsumoto, Ichiro Onoyama, Kazunori Imaizumi, Keiichi I. Nakayama

J Biol Chem. 2013 October 4; 288(40): 28488–28502. Published online 2013 August 16.
doi: 10.1074/jbc.M113.465179
PMCID: PMC3789950

59: Regulation of Stress-Inducible Phosphoprotein 1 Nuclear Retention by Protein Inhibitor of Activated STAT PIAS1

Iaci N. Soares, Fabiana A. Caetano, Jordan Pinder, Bruna Roz Rodrigues, Flavio H. Beraldo, Valeriy G. Ostapchenko, Chantal Durette, Grace Schenatto Pereira, Marilene H. Lopes, Nicolle Queiroz-Hazarbassanov, Isabela W. Cunha, Paulo I. Sanematsu, Sergio Suzuki, Luiz F. Bleggi-Torres, Caroline Schild-Poulter, Pierre Thibault, Graham Dellaire, Vilma R. Martins, Vania F. Prado, Marco A. M. Prado

Mol Cell Proteomics. 2013 November; 12(11): 3253–3270. Published online 2013 August 12.
doi: 10.1074/mcp.M113.031005
PMCID: PMC3820937

60: Human Mitochondrial Chaperone (mtHSP70) and Cysteine Desulfurase (NFS1) Bind Preferentially to the Disordered Conformation, Whereas Co-chaperone (HSC20) Binds to the Structured Conformation of the Iron-Sulfur Cluster Scaffold Protein (ISCU)

Kai Cai, Ronnie O. Frederick, Jin Hae Kim, Nichole M. Reinen, Marco Tonelli, John L. Markley
J Biol Chem. 2013 October 4; 288(40): 28755–28770. Published online 2013 August 12.
doi: 10.1074/jbc.M113.482042

PMCID: PMC3789972

61: Deubiquitinases Sharpen Substrate Discrimination during Membrane Protein Degradation from the ER

Zai-Rong Zhang, Juan S. Bonifacino, Ramanujan S. Hegde

Cell. 2013 August 1; 154(3): 609–622. doi: 10.1016/j.cell.2013.06.038
PMCID: PMC3732389

62: OTULIN restricts Met1-linked ubiquitination to control innate immune signaling

Berthe Katrine Fiil, Rune Busk Damgaard, Sebastian Alexander Wagner, Kirstin Keusekotten, Melanie Fritsch, Simon Bekker-Jensen, Niels Mailand, Chunaram Choudhary, David Komander, Mads Gyrd-Hansen

Mol Cell. Author manuscript; available in PMC 2014 October 13. Published in final edited form as: Mol Cell. 2013 June 27; 50(6): 818–830. doi: 10.1016/j.molcel.2013.06.004
PMCID: PMC4194427

63: Crystal structure of the Golgi casein kinase

Junyu Xiao, Vincent S. Tagliabracci, Jianzhong Wen, Soo-A Kim, Jack E. Dixon

Proc Natl Acad Sci U S A. 2013 June 25; 110(26): 10574–10579. Published online 2013 June 10. doi: 10.1073/pnas.1309211110

PMCID: PMC3696822

64: Mapping and Characterization of the Interaction Interface between Two Polypyrimidine-Tract Binding Proteins and a Nova-Type Protein of *Solanum tuberosum*

Shweta Shah, Nathaniel M. Butler, David J. Hannapel, A. Gururaj Rao

PLoS One. 2013; 8(5): e64783. Published online 2013 May 24.

doi: 10.1371/journal.pone.0064783

PMCID: PMC3663837

65: The Human CD8 β M-4 Isoform Dominant in Effector Memory T Cells Has Distinct Cytoplasmic Motifs That Confer Unique Properties

Deepshi Thakral, Maria M. Coman, Arunima Bandyopadhyay, Sunil Martin, James L. Riley, Paula B. Kavathas

PLoS One. 2013; 8(3): e59374. Published online 2013 March 22.

doi: 10.1371/journal.pone.0059374

PMCID: PMC3606432

66: Reversible inactivation of deubiquitinases by reactive oxygen species in vitro and in cells
Jin-Gu Lee, Kheewoong Baek, Nia Soetandyo, Yihong Ye

Nat Commun. 2013 March 5; 4: 1568. Published online 2013 March 5.

doi: 10.1038/ncomms2532

PMCID: PMC3615374

67: Evidence for an ABC-Type Riboflavin Transporter System in Pathogenic Spirochetes

Ranjit K. Deka, Chad A. Brautigam, Brent A. Bidy, Wei Z. Liu, Michael V. Norgard

mBio. 2013 Jan-Feb; 4(1): e00615-12. Published online 2013 February 12.

doi: 10.1128/mBio.00615-12

PMCID: PMC3573665

68: Proteasome inhibition prolongs survival during lethal hemorrhagic shock in rats

Harold H. Bach, IV, Heather M. LaPorte, Yee M. Wong, Richard L. Gamelli, Matthias Majetschak

J Trauma Acute Care Surg. Author manuscript; available in PMC 2014 October 2. Published in final edited form as: J Trauma Acute Care Surg. 2013 February; 74(2): 499–507.

doi: 10.1097/TA.0b013e31827d5db2

PMCID: PMC4182954

69: Refined Preparation and Use of Anti-diglycine Remnant (K- ϵ -GG) Antibody Enables Routine Quantification of 10,000s of Ubiquitination Sites in Single Proteomics Experiments

Namrata D. Udeshi, Tanya Svinkina, Philipp Mertins, Eric Kuhn, D. R. Mani, Jana W. Qiao, Steven A. Carr

Mol Cell Proteomics. 2013 March; 12(3): 825–831. Published online 2012 December 24.

doi: 10.1074/mcp.O112.027094

PMCID: PMC3591673

70: Transient kinetic analysis of USP2-catalyzed deubiquitination reveals a conformational rearrangement in the K48-linked diubiquitin substrate

William P. Bozza, Qin Liang, Ping Gong, Zhihao Zhuang

Biochemistry. Author manuscript; available in PMC 2013 December 18. Published in final edited form as: Biochemistry. 2012 December 18; 51(50): 10075–10086. Published online 2012

December 4. doi: 10.1021/bi3009104

PMCID: PMC3617563

71: Nonenzymatic Polymerization of Ubiquitin: Single-Step Synthesis and Isolation of Discrete Ubiquitin Oligomers

Vivian H. Trang, Ellen M. Valkevich, Shoko Minami, Yi-Chen Chen, Ying Ge, Eric R. Strieter
Angew Chem Int Ed Engl. Author manuscript; available in PMC 2014 July 7. Published in final edited form as: Angew Chem Int Ed Engl. 2012 December 21; 51(52): 13085–13088. Published online 2012 November 19. doi: 10.1002/anie.201207171

PMCID: PMC4083817

72: Involvement of WalK (VicK) Phosphatase Activity in Setting WalR (VicR) Response Regulator Phosphorylation Level and Limiting Crosstalk in *Streptococcus pneumoniae* D39 Cells

Kyle J. Wayne, Shuo Li, Krystyna M. Kazmierczak, Ho-Ching T. Tsui, Malcolm E. Winkler
Mol Microbiol. Author manuscript; available in PMC 2013 November 1. Published in final edited form as: Mol Microbiol. 2012 November; 86(3): 645–660. Published online 2012 September 27. doi: 10.1111/mmi.12006

PMCID: PMC3638944

73: TBK-1 promotes autophagy-mediated antimicrobial defense by controlling autophagosome maturation

Manohar Pilli, John Arko-Mensah, Marisa Ponpuak, Esteban Roberts, Sharon Master, Michael Mandell, Nicolas Dupont, Wojciech Ornatowski, Shanya Jiang, Steven Bradfute, Jack-Ansgar Bruun, Tom Egil Hansen, Terje Johansen, Vojo Deretic
Immunity. Author manuscript; available in PMC 2013 August 24. Published in final edited form as: Immunity. 2012 August 24; 37(2): 223–234. doi: 10.1016/j.immuni.2012.04.015

PMCID: PMC3428731

74: PDZ interactions regulate rapid turnover of the scaffolding protein EBP50 in microvilli
Damien Garbett, Anthony Bretscher

J Cell Biol. 2012 July 23; 198(2): 195–203. doi: 10.1083/jcb.201204008

PMCID: PMC3410424

75: The tyrosine phosphatase STEP: implications in schizophrenia and the molecular mechanism underlying antipsychotic medications

N C Carty, J Xu, P Kurup, J Brouillette, S M Goebel-Goody, D R Austin, P Yuan, G Chen, P R Correa, V Haroutunian, C Pittenger, P J Lombroso
Transl Psychiatry. 2012 July; 2(7): e137. Published online 2012 July 10.

doi: 10.1038/tp.2012.63

PMCID: PMC3410627

76: Molecular Basis of Lys-63-linked Polyubiquitination Inhibition by the Interaction between Human Deubiquitinating Enzyme OTUB1 and Ubiquitin-conjugating Enzyme UBC13

Yusuke Sato, Atsushi Yamagata, Sakurako Goto-Ito, Keiko Kubota, Rikako Miyamoto, Shinichiro Nakada, Shuya Fukai

J Biol Chem. 2012 July 27; 287(31): 25860–25868. Published online 2012 June 7.

doi: 10.1074/jbc.M112.364752

PMCID: PMC3406671

77: Characterization of a Glucosyltransferase Enzyme Involved in the Formation of Kaempferol and Quercetin Sphorosides in *Crocus sativus*

Almudena Trapero, Oussama Ahrazem, Angela Rubio-Moraga, Maria Luisa Jimeno, Maria Dolores Gómez, Lourdes Gómez-Gómez

Plant Physiol. 2012 August; 159(4): 1335–1354. Published online 2012 May 30.

doi: 10.1104/pp.112.198069

PMCID: PMC3425182

78: Monoubiquitination Promotes Calpain Cleavage of the Protein Phosphatase 2A (PP2A) Regulatory Subunit $\alpha 4$, Altering PP2A Stability and Microtubule-associated Protein Phosphorylation

Guy R. Watkins, Ning Wang, Matthew D. Mazalouskas, Rey J. Gomez, Chris R. Guthrie, Brian C. Kraemer, Susann Schweiger, Benjamin W. Spiller, Brian E. Wadzinski

J Biol Chem. 2012 July 13; 287(29): 24207–24215. Published online 2012 May 21.

doi: 10.1074/jbc.M112.368613

PMCID: PMC3397847

79: Role of Ubiquitin in Parainfluenza Virus 5 Particle Formation

Megan S. Harrison, Phuong Tieu Schmitt, Zifei Pei, Anthony P. Schmitt

J Virol. 2012 April; 86(7): 3474–3485. doi: 10.1128/JVI.06021-11

PMCID: PMC3302527

80: Toward Low-Cost Affinity Reagents: Lyophilized Yeast-scFv Probes Specific for Pathogen Antigens

Sean A. Gray, Kris M. Weigel, Ibne K. M. Ali, Annie A. Lakey, Jeremy Capalungan, Gonzalo J. Domingo, Gerard A. Cangelosi

PLoS One. 2012; 7(2): e32042. Published online 2012 February 20.

doi: 10.1371/journal.pone.0032042

PMCID: PMC3282784

81: Selectivity in Post-translational Biotin Addition to Five Human Carboxylases

Maria Ingaramo, Dorothy Beckett

J Biol Chem. 2012 January 13; 287(3): 1813–1822. Published online 2011 November 28.

doi: 10.1074/jbc.M111.275982

PMCID: PMC3265863

82: A mechanism for tunable autoinhibition in the structure of a human Ca^{2+} /calmodulin-dependent kinase II holoenzyme

Luke H. Chao, Margaret M. Stratton, Il-Hyung Lee, Oren S. Rosenberg, Joshua Levitz, Daniel J. Mandell, Tanja Kortemme, Jay T. Groves, Howard Schulman, John Kuriyan

Cell. Author manuscript; available in PMC 2012 March 2. Published in final edited form as: Cell. 2011 September 2; 146(5): 732–745. doi: 10.1016/j.cell.2011.07.038 Correction in: volume 147 on page 704
PMCID: PMC3184253

83: Eliminating SF-1 (NR5A1) Sumoylation In Vivo Results in Ectopic Hedgehog Signaling and Disruption of Endocrine Development
Florence Y. Lee, Emily J. Faivre, Miyuki Suzawa, Erik Lontok, Daniel Ebert, Fang Cai, Denise D. Belsham, Holly A. Ingraham
Dev Cell. Author manuscript; available in PMC 2012 August 16. Published in final edited form as: Dev Cell. 2011 August 16; 21(2): 315–327. Published online 2011 August 4.
doi: 10.1016/j.devcel.2011.06.028
PMCID: PMC3157481

84: Immunity-related GTPase M (IRGM) Proteins Influence the Localization of Guanylate-binding Protein 2 (GBP2) by Modulating Macroautophagy
Maria K. Traver, Stanley C. Henry, Viviana Cantillana, Tim Oliver, Julia P. Hunn, Jonathan C. Howard, Sandra Beer, Klaus Pfeffer, Jörn Coers, Gregory A. Taylor
J Biol Chem. 2011 September 2; 286(35): 30471–30480. Published online 2011 July 12.
doi: 10.1074/jbc.M111.251967 Correction in: J Biol Chem. 2013 April 19; 288(16): 11504.
PMCID: PMC3162407

85: The variable detergent sensitivity of proteases that are utilized for recombinant protein affinity tag removal
James M. Vergis, Michael C. Wiener
Protein Expr Purif. Author manuscript; available in PMC 2012 August 1. Published in final edited form as: Protein Expr Purif. 2011 August; 78(2): 139–142. Published online 2011 April 24. doi: 10.1016/j.pep.2011.04.011
PMCID: PMC3130611

86: 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, Benjamin W. Spiller
J Biol Chem. 2011 May 20; 286(20): 17665–17671. Published online 2011 March 29.
doi: 10.1074/jbc.M111.222414
PMCID: PMC3093842

87: Biotinylation, a Post-translational Modification Controlled by the Rate of Protein-Protein Association
Maria Ingaramo, Dorothy Beckett
J Biol Chem. 2011 April 15; 286(15): 13071–13078. Published online 2011 February 22.
doi: 10.1074/jbc.M110.183624
PMCID: PMC3075653

88: Challenges associated with the targeted delivery of gelonin to claudin-expressing cancer cells with the use of activatable cell penetrating peptides to enhance potency

Xiaoqin Yuan, Xinjian Lin, Gerald Manorek, Stephen B Howell

BMC Cancer. 2011; 11: 61. Published online 2011 February 8. doi: 10.1186/1471-2407-11-61
PMCID: PMC3042419

89: SUMOylation of DLX3 by SUMO1 promotes its transcriptional activity

O. Duverger, S.X. Chen, D. Lee, T. Li, P.B. Chock, M.I. Morasso

J Cell Biochem. Author manuscript; available in PMC 2012 February 1. Published in final edited form as: J Cell Biochem. 2011 February; 112(2): 445–452. doi: 10.1002/jcb.22891
PMCID: PMC3180851

90: A Quantitative Model of Thermal Stabilization and Destabilization of Proteins by Ligands
Piotras Cimmperman, Lina Baranauskienė, Simona Jachimovičiūtė, Jelena Jachno, Jolanta Torresan, Vilma Michailovienė, Jurgita Matulienė, Jolanta Sereikaitė, Vladas Bumelis, Daumantas Matulis

Biophys J. 2008 October 1; 95(7): 3222–3231. Published online 2008 July 3.
doi: 10.1529/biophysj.108.134973
PMCID: PMC2547457

91: Characterization of ubiquitin and ubiquitin-like-protein isopeptidase activities

Benjamin Nicholson, Craig A. Leach, Seth J. Goldenberg, Dana M. Francis, Matthew P.

Kodrasov, Xufan Tian, John Shanks, David E. Sterner, Alejandro Bernal, Michael R. Mattern, Keith D. Wilkinson, Tauseef R. Butt

Protein Sci. 2008 June; 17(6): 1035–1043. doi: 10.1110/ps.083450408
PMCID: PMC2386736

92: SENP1-mediated GATA1 deSUMOylation is critical for definitive erythropoiesis

Luyang Yu, Weidong Ji, Haifeng Zhang, Matthew J. Renda, Yun He, Sharon Lin, Ee-chun Cheng, Hong Chen, Diane S. Krause, Wang Min

J Exp Med. 2010 June 7; 207(6): 1183–1195. doi: 10.1084/jem.20092215
PMCID: PMC2882842

93: Crystallization and preliminary crystallographic studies of the catalytic subunits of human pyruvate dehydrogenase phosphatase isoforms 1 and 2

Junko Kato, Masato Kato

Acta Crystallogr Sect F Struct Biol Cryst Commun. 2010 March 1; 66(Pt 3): 342–345.

Published online 2010 February 27. doi: 10.1107/S1744309110003131
PMCID: PMC2833053

94: Kinetic Characterization of the WalRKSpn (VicRK) Two-Component System of *Streptococcus pneumoniae*: Dependence of WalKSpn (VicK) Phosphatase Activity on Its PAS Domain

Alina D. Gutu, Kyle J. Wayne, Lok-To Sham, Malcolm E. Winkler

J Bacteriol. 2010 May; 192(9): 2346–2358. Published online 2010 February 26.
doi: 10.1128/JB.01690-09

PMCID: PMC2863487

95: Identification of SRC3/AIB1 as a Preferred Coactivator for Hormone-activated Androgen Receptor

X. Edward Zhou, Kelly M. Suino-Powell, Jun Li, Yuanzheng He, Jeffrey P. MacKeigan, Karsten Melcher, Eu-Leong Yong, H. Eric Xu

J Biol Chem. 2010 March 19; 285(12): 9161–9171. Published online 2010 January 19.

doi: 10.1074/jbc.M109.085779

PMCID: PMC2838335

96: Cerebrospinal Fluid Secretory Ca²⁺-Dependent Phospholipase A2 Activity Is Increased in Alzheimer Disease

Sonia Chalbot, Henrik Zetterberg, Kaj Blennow, Tormod Fladby, Inge Grundke-Iqbal, Khalid Iqbal

Clin Chem. Author manuscript; available in PMC 2010 April 21. Published in final edited form as: Clin Chem. 2009 December; 55(12): 2171–2179. Published online 2009 October 22.

doi: 10.1373/clinchem.2009.130286

PMCID: PMC2857971

97: The molecular basis for the regulation of the CBC by the importins

Sandra M.G. Dias, Kristin F. Wilson, Katherine S. Rojas, Andre L.B. Ambrosio, Richard A. Cerione

Nat Struct Mol Biol. Author manuscript; available in PMC 2010 March 1. Published in final edited form as: Nat Struct Mol Biol. 2009 September; 16(9): 930–937. Published online 2009 August 9. doi: 10.1038/nsmb.1649

PMCID: PMC2782468

98: Molecular Identification and Expression Analysis of Filaggrin-2, a Member of the S100 Fused-Type Protein Family

Zhihong Wu, Britta Hansmann, Ulf Meyer-Hoffert, Regine Gläser, Jens-Michael Schröder

PLoS ONE. 2009; 4(4): e5227. Published online 2009 April 22.

doi: 10.1371/journal.pone.0005227

PMCID: PMC2668185

99: Identification of Lympho-Epithelial Kazal-Type Inhibitor 2 in Human Skin as a Kallikrein-Related Peptidase 5-Specific Protease Inhibitor

Ulf Meyer-Hoffert, Zhihong Wu, Jens-Michael Schröder

PLoS ONE. 2009; 4(2): e4372. Published online 2009 February 3.

doi: 10.1371/journal.pone.0004372

PMCID: PMC2631147

100: The crystal structure of the catalytic domain of a eukaryotic guanylate cyclase

Jonathan A Winger, Emily R Derbyshire, Meindert H Lamers, Michael A Marletta, John Kuriyan

BMC Struct Biol. 2008; 8: 42. Published online 2008 October 7. doi: 10.1186/1472-6807-8-42

PMCID: PMC2576301

101: Decreased Recognition of SUMO-Sensitive Target Genes following Modification of SF-1 (NR5A1)

Lioudmila A. Campbell, Emily J. Faivre, Matthew D. Show, Jared G. Ingraham, Jeremy Flinders, John D. Gross, Holly A. Ingraham

Mol Cell Biol. 2008 December; 28(24): 7476–7486. Published online 2008 October 6.

doi: 10.1128/MCB.00103-08

PMCID: PMC2593425

102: Enhanced protein expression in the baculovirus/insect cell system using engineered SUMO fusions

Li Liu, Joshua Spurrier, Tauseef R. Butt, James E. Strickler

Protein Expr Purif. Author manuscript; available in PMC 2009 November 1. Published in final edited form as: Protein Expr Purif. 2008 November; 62(1): 21–28. Published online 2008 August 5. doi: 10.1016/j.pep.2008.07.010

PMCID: PMC2585507

103: Nucleotidylation of the VPg Protein of a Human Norovirus by its Proteinase-Polymerase Precursor Protein

Gaël Belliot, Stanislav V. Sosnovtsev, Kyeong-Ok Chang, Peter McPhie, Kim Y. Green

Virology. Author manuscript; available in PMC 2009 April 25. Published in final edited form as: Virology. 2008 April 25; 374(1): 33–49. Published online 2008 January 30.

doi: 10.1016/j.virol.2007.12.028

PMCID: PMC2386983

104: Structural transition in Bcl-xL and its potential association with mitochondrial calcium ion transport

Sreekanth Rajan, Minjoo Choi, Quoc Toan Nguyen, Hong Ye, Wei Liu, Hui Ting Toh, CongBao Kang, Neelagandan Kamariah, Chi Li, Huiya Huang, Carl White, Kwanghee Baek, Gerhard Grüber, Ho Sup Yoon

Sci Rep. 2015; 5: 10609. Published online 2015 May 29. doi: 10.1038/srep10609

PMCID: PMC4448555

105: K63 linked ubiquitin chain formation is a signal for HIF1A degradation by Chaperone-Mediated Autophagy

Joao Vasco Ferreira, Ana Rosa Soares, Jose Silva Ramalho, Paulo Pereira, Henrique Girao

Sci Rep. 2015; 5: 10210. Published online 2015 May 11. doi: 10.1038/srep10210

PMCID: PMC4426689

106: Oligomerization of FtsZ converts the FtsZ tail motif (CCTP) into a multivalent ligand with high avidity for partners ZipA and SImA

Shishen Du, Kyung-Tae Park, Joe Lutkenhaus

Mol Microbiol. Author manuscript; available in PMC 2015 March 18. Published in final edited form as: Mol Microbiol. 2015 January; 95(2): 173–188. Published online 2014 December 19.

doi: 10.1111/mmi.12854

PMCID: PMC4364296

107: Catalytic Function and Substrate Specificity of the Papain-Like Protease Domain of nsp3 from the Middle East Respiratory Syndrome Coronavirus

Yahira M. Báez-Santos, Anna M. Mielech, Xufang Deng, Susan Baker, Andrew D. Mesecar
J Virol. 2014 November; 88(21): 12511–12527. doi: 10.1128/JVI.01294-14
 PMID: PMC4248884

108: Recruitment of Cbl-b to B Cell Antigen Receptor Couples Antigen Recognition to Toll-Like Receptor 9 Activation in Late Endosomes

Margaret Veselits, Azusa Tanaka, Stanley Lipkowitz, Shannon O'Neill, Roger Sciammas, Alison Finnegan, Jian Zhang, Marcus R. Clark
PLoS One. 2014; 9(3): e89792. Published online 2014 March 20.
 doi: 10.1371/journal.pone.0089792
 PMID: PMC3961229

109: 19F NMR Reveals Multiple Conformations at the Dimer Interface of the Non-Structural Protein 1 Effector Domain from Influenza A Virus

James M. Aramini, Keith Hamilton, Li-Chung Ma, G. V. T. Swapna, Paul G. Leonard, John E. Ladbury, Robert M. Krug, Gaetano T. Montelione
Structure. Author manuscript; available in PMC 2015 April 8. Published in final edited form as: *Structure.* 2014 April 8; 22(4): 515–525. Published online 2014 February 27.
 doi: 10.1016/j.str.2014.01.010
 PMID: PMC4110948

110: Induction of Gsk3 β - β -TrCP Interaction Is Required for Late Phase Stabilization of β -Catenin in Canonical Wnt Signaling

Chenxi Gao, Guangming Chen, Guillermo Romero, Stergios Moschos, Xiang Xu, Jing Hu
J Biol Chem. 2014 March 7; 289(10): 7099–7108. Published online 2014 January 22.
 doi: 10.1074/jbc.M113.532606
 PMID: PMC3945370

111: Ubiquitination-dependent quality control of hERG K⁺ channel with acquired and inherited conformational defect at the plasma membrane

Pirjo M. Apaja, Brian Foo, Tsukasa Okiyoneda, William C. Valinsky, Herve Barriere, Roxana Atanasiu, Eckhard Ficker, Gergely L. Lukacs, Alvin Shrier
Mol Biol Cell. 2013 December 15; 24(24): 3787–3804. doi: 10.1091/mbc.E13-07-0417
 PMID: PMC3861077

112: The tails of apical scaffolding proteins EBP50 and E3KARP regulate their localization and dynamics

Damien Garbett, Cécile Sauvanet, Raghuvir Viswanatha, Anthony Bretscher
Mol Biol Cell. 2013 November 1; 24(21): 3381–3392. doi: 10.1091/mbc.E13-06-0330
 PMID: PMC3814156

113: Error rates for nanopore discrimination among cytosine, methylcytosine, and hydroxymethylcytosine along individual DNA strands

Jacob Schreiber, Zachary L. Wescoe, Robin Abu-Shumays, John T. Vivian, Baldandorj Baatar, Kevin Karplus, Mark Akeson

Proc Natl Acad Sci U S A. 2013 November 19; 110(47): 18910–18915. Published online 2013 October 28. doi: 10.1073/pnas.1310615110

PMCID: PMC3839712

114: Structural and Functional Analyses of DNA-Sensing and Immune Activation by Human cGAS

Kazuki Kato, Ryohei Ishii, Eiji Goto, Ryuichiro Ishitani, Fuminori Tokunaga, Osamu Nureki
PLoS One. 2013; 8(10): e76983. Published online 2013 October 7.

doi: 10.1371/journal.pone.0076983

PMCID: PMC3792152

115: Antimicrobial Activity of High-Mobility-Group Box 2: a New Function to a Well-Known Protein

Robert K uchler, Bjoern O. Schroeder, Simon U. Jaeger, Eduard F. Stange, Jan Wehkamp

Antimicrob Agents Chemother. 2013 October; 57(10): 4782–4793. doi: 10.1128/AAC.00805-13

PMCID: PMC3811396

116: Using Antiubiquitin Antibodies to Probe the Ubiquitination State Within rhTRIM5 α Cytoplasmic Bodies

Cindy M. Danielson, Thomas J. Hope

AIDS Res Hum Retroviruses. 2013 October; 29(10): 1373–1385. doi: 10.1089/aid.2013.0029

PMCID: PMC3785812

117: Jun Activation Domain-binding Protein 1 (JAB1) Is Required for the Optimal Response to Interferons

Ryuta Muromoto, Maiko Nakajima, Koki Hirashima, Toru Hirao, Shigeyuki Kon, Kazuya Shimoda, Kenji Oritani, Tadashi Matsuda

J Biol Chem. 2013 October 25; 288(43): 30969–30979. Published online 2013 September 16.

doi: 10.1074/jbc.M113.485847

PMCID: PMC3829410

118: SARM1 and TRAF6 bind to and stabilize PINK1 on depolarized mitochondria

Hitoshi Murata, Masakiyo Sakaguchi, Ken Kataoka, Nam-ho Huh

Mol Biol Cell. 2013 September 15; 24(18): 2772–2784. doi: 10.1091/mbc.E13-01-0016

PMCID: PMC3771941

119: Measuring activity in the ubiquitin-proteasome system: From large scale discoveries to single cells analysis

Adam T. Melvin, Gregory S. Woss, Jessica H. Park, Marcey L. Waters, Nancy L. Allbritton

Cell Biochem Biophys. Author manuscript; available in PMC 2014 September 1. Published in final edited form as: Cell Biochem Biophys. 2013 September; 67(1): 75–89.

doi: 10.1007/s12013-013-9621-9

PMCID: PMC3758771

120: Redox and Light Control the Heme-Sensing Activity of AppA

Liang Yin, Vladimira Dragnea, George Feldman, Loubna A. Hammad, Jonathan A. Karty, Charles E. Dann, III, Carl E. Bauer

mBio. 2013 Sep-Oct; 4(5): e00563-13. Published online 2013 August 27.

doi: 10.1128/mBio.00563-13

PMCID: PMC3760249

121: Stabilization of the Transcription Factor Foxp3 by the Deubiquitinase USP7 Increases Treg-Cell-Suppressive Capacity

Jorg van Loosdregt, Veerle Fleskens, Juan Fu, Arjan B. Brenkman, Cornelis P.J. Bekker, Cornelie E.G.M. Pals, Jenny Meerding, Celia R. Berkers, Joseph Barbi, Andrea Gröne, Alice J.A.M. Sijts, Madelon M. Maurice, Eric Kalkhoven, Berent J. Prakken, Huib Ovaa, Fan Pan, Dietmar M.W. Zaiss, Paul J. Coffey

Immunity. Author manuscript; available in PMC 2014 August 15. Published in final edited form as: Immunity. 2013 August 22; 39(2): 259–271. doi: 10.1016/j.immuni.2013.05.018

PMCID: PMC4133784

122: E2 ubiquitin conjugating enzymes regulate the deubiquitinating activity of OTUB1

Reuven Wiener, Anthony T. DiBello, Patrick Lombardi, Catherine M. Guzzo, Xiangbin Zhang, Michael J. Matunis, Cynthia Wolberger

Nat Struct Mol Biol. Author manuscript; available in PMC 2014 March 4. Published in final edited form as: Nat Struct Mol Biol. 2013 September; 20(9): 1033–1039. Published online 2013 August 18. doi: 10.1038/nsmb.2655

PMCID: PMC3941643

123: Haxe1 controls ROS generation of vertebrate Rac1-dependent NADPH oxidase complexes

Mads Daugaard, Roberto Nitsch, Babak Razaghi, Lindsay McDonald, Ameer Jarrar, Stéphanie Torino, Sonia Castillo-Lluva, Barak Rotblat, Liheng Li, Angeliki Malliri, Emmanuel Lemichez, Amel Mettouchi, Jason N. Berman, Josef M. Penninger, Poul H. Sorensen

Nat Commun. 2013 July 17; 4: 2180. Published online 2013 July 17. doi: 10.1038/ncomms3180

PMCID: PMC3759041

124: Human Parainfluenza Virus Type 2 V Protein Inhibits TRAF6-Mediated Ubiquitination of IRF7 To Prevent TLR7- and TLR9-Dependent Interferon Induction

Yoshinori Kitagawa, Mayu Yamaguchi, Min Zhou, Machiko Nishio, Masae Itoh, Bin Gotoh

J Virol. 2013 July; 87(14): 7966–7976. doi: 10.1128/JVI.03525-12

PMCID: PMC3700191

125: Pressure-Induced Endocytic Degradation of the Saccharomyces cerevisiae Low-Affinity Tryptophan Permease Tat1 Is Mediated by Rsp5 Ubiquitin Ligase and Functionally Redundant PPxY Motif Proteins

Asaha Suzuki, Takahiro Mochizuki, Satoshi Uemura, Toshiki Hiraki, Fumiyoshi Abe

Eukaryot Cell. 2013 July; 12(7): 990–997. doi: 10.1128/EC.00049-13

PMCID: PMC3697464

126: MOR Is Not Enough: Identification of Novel mu-Opioid Receptor Interacting Proteins Using Traditional and Modified Membrane Yeast Two-Hybrid Screens
Jessica Petko, Stephanie Justice-Bitner, Jay Jin, Victoria Wong, Saranya Kittanakom, Thomas N. Ferraro, Igor Stagljjar, Robert Levenson
PLoS One. 2013; 8(6): e67608. Published online 2013 June 28.
doi: 10.1371/journal.pone.0067608
PMCID: PMC3695902

127: Decoupling the Role of Ubiquitination for the Dislocation Versus Degradation of Major Histocompatibility Complex (MHC) Class I Proteins during Endoplasmic Reticulum-associated Degradation (ERAD)
Xiaoli Wang, Y. Y. Lawrence Yu, Nancy Myers, Ted H. Hansen
J Biol Chem. 2013 August 9; 288(32): 23295–23306. Published online 2013 June 25.
doi: 10.1074/jbc.M113.482018
PMCID: PMC3743500

128: 53BP1 is a reader of the DNA damage-induced H2A Lys15 ubiquitin mark
Amélie Fradet-Turcotte, Marella D. Canny, Cristina Escribano-Díaz, Alexandre Orthwein, Charles C.Y. Leung, Hao Huang, Marie-Claude Landry, Julianne Kitevski-LeBlanc, Sylvie M. Noordermeer, Frank Sicheri, Daniel Durocher
Nature. Author manuscript; available in PMC 2014 March 15. Published in final edited form as: Nature. 2013 July 4; 499(7456): 50–54. Published online 2013 June 12.
doi: 10.1038/nature12318
PMCID: PMC3955401

129: Stabilization of the μ -Opioid Receptor by Truncated Single Transmembrane Splice Variants through a Chaperone-like Action
Jin Xu, Ming Xu, Taylor Brown, Grace C. Rossi, Yasmin L. Hurd, Charles E. Inturrisi, Gavril W. Pasternak, Ying-Xian Pan
J Biol Chem. 2013 July 19; 288(29): 21211–21227. Published online 2013 June 11.
doi: 10.1074/jbc.M113.458687
PMCID: PMC3774389

130: A Novel Role of the N Terminus of Budding Yeast Histone H3 Variant Cse4 in Ubiquitin-Mediated Proteolysis
Wei Chun Au, Anthony R. Dawson, David W. Rawson, Sara B. Taylor, Richard E. Baker, Munira A. Basrai
Genetics. 2013 June; 194(2): 513–518. doi: 10.1534/genetics.113.149898
PMCID: PMC3664860

131: NEW TOOLS FOR THE BRAINBOW TOOLBOX
Dawen Cai, Kimberly B. Cohen, Tuanlian Luo, Jeff W. Lichtman, Joshua R. Sanes
Nat Methods. Author manuscript; available in PMC 2013 November 5. Published in final edited form as: Nat Methods. 2013 May 5; 10(6): 540–547. Published online 2013 May 5.
doi: 10.1038/nmeth.2450
PMCID: PMC3713494

132: Preliminary X-ray crystallographic studies of an N-terminal domain of unknown function from a putative glycosyltransferase from *Streptococcus parasanguinis*

Hua Zhang, Fan Zhu, Lei Ding, Meixian Zhou, Ren Wu, Hui Wu

Acta Crystallogr Sect F Struct Biol Cryst Commun. 2013 May 1; 69(Pt 5): 520–523. Published online 2013 April 30. doi: 10.1107/S1744309113007161

PMCID: PMC3660891

133: Analysis of SUMOylated proteins using SUMO-traps

Elisa Da Silva-Ferrada, Wendy Xolalpa, Valérie Lang, Fabienne Aillet, Itziar Martin-Ruiz, Carlos F. de la Cruz-Herrera, Fernando Lopitz-Otsoa, Arkaitz Carracedo, Seth J. Goldenberg, Carmen Rivas, Patrick England, Manuel S. Rodríguez

Sci Rep. 2013; 3: 1690. Published online 2013 April 22. doi: 10.1038/srep01690

PMCID: PMC3631770

134: Single-molecule sorting reveals how ubiquitylation affects substrate recognition and activities of FBH1 helicase

Tokiha Masuda-Ozawa, Trish Hoang, Yeon-Soo Seo, Lin-Feng Chen, Maria Spies

Nucleic Acids Res. 2013 April; 41(6): 3576–3587. Published online 2013 February 6.

doi: 10.1093/nar/gkt056

PMCID: PMC3616717

135: Crystal structures of two phytohormone signal-transducing α/β hydrolases: karrikin-signaling KAI2 and strigolactone-signaling DWARF14

Li-Hua Zhao, X Edward Zhou, Zhong-Shan Wu, Wei Yi, Yong Xu, Suling Li, Ting-Hai Xu, Yue Liu, Run-Ze Chen, Amanda Kovach, Yangyong Kang, Li Hou, Yuanzheng He, Cen Xie, Wanling Song, Dafang Zhong, Yechun Xu, Yonghong Wang, Jiayang Li, Chenghai Zhang, Karsten Melcher, H Eric Xu

Cell Res. 2013 March; 23(3): 436–439. Published online 2013 February 5.

doi: 10.1038/cr.2013.19

PMCID: PMC3587710

136: Unfoldase-mediated protein translocation through an α -hemolysin nanopore

Jeff Nivala, Douglas B Marks, Mark Akeson

Nat Biotechnol. Author manuscript; available in PMC 2013 September 13. Published in final edited form as: Nat Biotechnol. 2013 March; 31(3): 247–250. Published online 2013 February 3.

doi: 10.1038/nbt.2503

PMCID: PMC3772521

137: Proteasomal degradation of herpes simplex virus capsids in macrophages releases DNA to the cytosol for recognition by DNA sensors

Kristy A. Horan, Kathrine Hansen, Martin R. Jakobsen, Christian K. Holm, Stine Søby, Leonie Unterholzner, Mikayla Thompson, John A. West, Marie B. Iversen, Simon B. Rasmussen, Svend Ellermann-Eriksen, Evelyn Kurt-Jones, Santo Landolfo, Blossom Damania, Jesper Melchjorsen, Andrew G. Bowie, Katherine A. Fitzgerald, Søren R. Paludan

J Immunol. Author manuscript; available in PMC 2014 March 1. Published in final edited form as: J Immunol. 2013 March 1; 190(5): 2311–2319. Published online 2013 January 23.

doi: 10.4049/jimmunol.1202749

PMCID: PMC3578088

138: Trypanosoma brucei S-Adenosylmethionine Decarboxylase N Terminus Is Essential for Allosteric Activation by the Regulatory Subunit Prozyme

Nahir Velez, Chad A. Brautigam, Margaret A. Phillips

J Biol Chem. 2013 February 15; 288(7): 5232–5240. Published online 2013 January 3.

doi: 10.1074/jbc.M112.442475

PMCID: PMC3576127

139: MG53-induced IRS-1 ubiquitination negatively regulates skeletal myogenesis and insulin signaling

Jae-Sung Yi, Jun Sub Park, Young-Mi Ham, Nga Nguyen, Na-Rae Lee, Jin Hong, Bong-Woo Kim, Hyun Lee, Chang-Seok Lee, Byung-Cheon Jeong, Hyun Kyu Song, Hana Cho, Yoon Ki Kim, Jae-Seon Lee, Kyong-Soo Park, Hakseob Shin, Inho Choi, Seung Hee Lee, Woo Jin Park, Shi-Young Park, Cheol Soo Choi, Peihui Lin, Malith Karunasiri, Tao Tan, Pu Duann, Hua Zhu, Jianjie Ma, Young-Gyu Ko

Nat Commun. Author manuscript; available in PMC 2014 March 4. Published in final edited form as: Nat Commun. 2013; 4: 2354. doi: 10.1038/ncomms3354

PMCID: PMC3941707

140: Small ubiquitin-like modifying protein isopeptidase assay based on poliovirus RNA polymerase activity

Jamie J. Arnold, Alejandro Bernal, Uzo Uche, David E. Sterner, Tauseef R. Butt, Craig E. Cameron, Michael R. Mattern

Anal Biochem. Author manuscript; available in PMC 2007 November 26. Published in final edited form as: Anal Biochem. 2006 March 15; 350(2): 214–221. Published online 2005 November 17. doi: 10.1016/j.ab.2005.11.001

PMCID: PMC2094218

141: Androgen receptor degradation by the E3 ligase CHIP modulates mitotic arrest in prostate cancer cells

S Sarkar, DL Brautigam, SJ Parsons, JM Larner

Oncogene. Author manuscript; available in PMC 2014 January 23. Published in final edited form as: Oncogene. 2014 January 2; 33(1): 26–33. Published online 2012 December 17.

doi: 10.1038/onc.2012.561

PMCID: PMC3899837

142: The Interaction of Prp2 with a Defined Region of the Intron Is Required for the First Splicing Reaction

Hsueh-Lien Liu, Soo-Chen Cheng

Mol Cell Biol. 2012 December; 32(24): 5056–5066. doi: 10.1128/MCB.01109-12

PMCID: PMC3510548

143: The Ω -loop lid domain of phosphoenolpyruvate carboxykinase is essential for catalytic function

Troy A. Johnson, Todd Holyoak

Biochemistry. Author manuscript; available in PMC 2013 November 27. Published in final edited form as: Biochemistry. 2012 November 27; 51(47): 9547–9559. Published online 2012

November 14. doi: 10.1021/bi301278t

PMCID: PMC3525453

144: MultiDsk: A Ubiquitin-Specific Affinity Resin

Marcus D. Wilson, Marco Saponaro, Mathias A. Leidl, Jesper Q. Svejstrup

PLoS One. 2012; 7(10): e46398. Published online 2012 October 3.

doi: 10.1371/journal.pone.0046398

PMCID: PMC3463603

145: Non-transcriptional Priming and Deubiquitination Regulate NLRP3 Inflammasome Activation

Christine Juliana, Teresa Fernandes-Alnemri, Seokwon Kang, Andrew Farias, Fengsong Qin, Emad S. Alnemri

J Biol Chem. 2012 October 19; 287(43): 36617–36622. Published online 2012 September 4.

doi: 10.1074/jbc.M112.407130

PMCID: PMC3476327

146: Polyubiquitin Is Required for Growth, Development and Pathogenicity in the Rice Blast Fungus *Magnaporthe oryzae*

Yeonyee Oh, William L. Franck, Sang-Oh Han, Angela Shows, Emine Gokce, David C.

Muddiman, Ralph A. Dean

PLoS One. 2012; 7(8): e42868. Published online 2012 August 10.

doi: 10.1371/journal.pone.0042868

PMCID: PMC3416782

147: Cathepsin G-regulated Release of Formyl Peptide Receptor Agonists Modulate Neutrophil Effector Functions

Josh C. Woloszynek, Ying Hu, Christine T. N. Pham

J Biol Chem. 2012 October 5; 287(41): 34101–34109. Published online 2012 August 9.

doi: 10.1074/jbc.M112.394452

PMCID: PMC3464519

148: Activity of the tetrapyrrole regulator CrtJ is controlled by oxidation of a redox active cysteine located in the DNA binding domain

Zhuo Cheng, Jiang Wu, Aaron Setterdahl, Khalilah Reddie, Kate Carroll, Loubna A. Hammad, Jonathan A. Karty, Carl E. Bauer

Mol Microbiol. Author manuscript; available in PMC 2013 August 1. Published in final edited form as: Mol Microbiol. 2012 August; 85(4): 734–746. Published online 2012 July 16.

doi: 10.1111/j.1365-2958.2012.08135.x

PMCID: PMC3418406

149: TP-RT Domain Interactions of Duck Hepatitis B Virus Reverse Transcriptase in cis and in trans during Protein-Primed Initiation of DNA Synthesis In Vitro

Rajeev K. Boregowda, Christina Adams, Jianming Hu

J Virol. 2012 June; 86(12): 6522–6536. doi: 10.1128/JVI.00086-12

PMCID: PMC3393569

150: Functional Dissection of N-Acetylglutamate Synthase (ArgA) of *Pseudomonas aeruginosa* and Restoration of Its Ancestral N-Acetylglutamate Kinase Activity

Enea Sancho-Vaello, María L. Fernández-Murga, Vicente Rubio

J Bacteriol. 2012 June; 194(11): 2791–2801. doi: 10.1128/JB.00125-12

PMCID: PMC3370605

151: Specific Residues of a Conserved Domain in the N Terminus of the Human Cytomegalovirus pUL50 Protein Determine Its Intranuclear Interaction with pUL53

Jens Milbradt, Sabrina Auerochs, Madhumati Sevana, Yves A. Muller, Heinrich Sticht, Manfred Marschall

J Biol Chem. 2012 July 6; 287(28): 24004–24016. Published online 2012 May 15.

doi: 10.1074/jbc.M111.331207

PMCID: PMC3390675

152: In Vitro Epsilon RNA-Dependent Protein Priming Activity of Human Hepatitis B Virus Polymerase

Scott A. Jones, Rajeev Boregowda, Thomas E. Spratt, Jianming Hu

J Virol. 2012 May; 86(9): 5134–5150. doi: 10.1128/JVI.07137-11 Correction in: J Virol. 2013 November; 87(22): 12504.

PMCID: PMC3347376

153: Structural and Biochemical Characterization of Human Mitochondrial Branched-chain α -Ketoacid Dehydrogenase Phosphatase

R. Max Wynn, Jun Li, Chad A. Brautigam, Jacinta L. Chuang, David T. Chuang

J Biol Chem. 2012 March 16; 287(12): 9178–9192. Published online 2012 January 30.

doi: 10.1074/jbc.M111.314963

PMCID: PMC3308798

154: Analysis of oxygen/glucose-deprivation-induced changes in SUMO3 conjugation using SILAC-based quantitative proteomics

Wei Yang, J. Will Thompson, Zhengfeng Wang, Liangli Wang, Huaxin Sheng, Matthew W. Foster, M. Arthur Moseley, Wulf Paschen

J Proteome Res. Author manuscript; available in PMC 2013 April 17. Published in final edited form as: J Proteome Res. 2012 February 3; 11(2): 1108–1117. Published online 2011 December 1. doi: 10.1021/pr200834f

PMCID: PMC3628696

155: Ubiquitin Ligase Substrate Identification through Quantitative Proteomics at Both the Protein and Peptide Levels

Kimberly A. Lee, Lisa P. Hammerle, Paul S. Andrews, Matthew P. Stokes, Tomas Mustelin, Jeffrey C. Silva, Roy A. Black, John R. Doedens
J Biol Chem. 2011 December 2; 286(48): 41530–41538. Published online 2011 October 10.
doi: 10.1074/jbc.M111.248856
PMCID: PMC3308864

156: The MHC I immunopeptidome conveys to the cell surface an integrative view of cellular regulation
Etienne Caron, Krystel Vincent, Marie-Hélène Fortier, Jean-Philippe Laverdure, Alexandre Bramoullé, Marie-Pierre Hardy, Grégory Voisin, Philippe P Roux, Sébastien Lemieux, Pierre Thibault, Claude Perreault
Mol Syst Biol. 2011; 7: 533. Published online 2011 September 27. doi: 10.1038/msb.2011.68
PMCID: PMC3202804

157: Crystal Structures of Histone and p53 Methyltransferase SmyD2 Reveal a Conformational Flexibility of the Autoinhibitory C-Terminal Domain
Yuanyuan Jiang, Nualpun Sirinupong, Joseph Brunzelle, Zhe Yang
PLoS One. 2011; 6(6): e21640. Published online 2011 June 28.
doi: 10.1371/journal.pone.0021640
PMCID: PMC3125274

158: The Ubiquitin-Specific Protease USP34 Regulates Axin Stability and Wnt/ β -Catenin Signaling
Tony T. H. Lui, Celine Lacroix, Syed M. Ahmed, Seth J. Goldenberg, Craig A. Leach, Avais M. Daulat, Stephane Angers
Mol Cell Biol. 2011 May; 31(10): 2053–2065. doi: 10.1128/MCB.01094-10
PMCID: PMC3133363

159: Comparison of SUMO fusion technology with traditional gene fusion systems: Enhanced expression and solubility with SUMO
Jeffrey G. Marblestone, Suzanne C. Edavettal, Yiting Lim, Peter Lim, Xun Zuo, Tauseef R. Butt
Protein Sci. 2006 January; 15(1): 182–189. doi: 10.1110/ps.051812706
PMCID: PMC2242369

160: Immunization with a ZmpB-Based Protein Vaccine Could Protect against Pneumococcal Diseases in Mice
Yi Gong, Wenchun Xu, Yali Cui, Xuemei Zhang, Run Yao, Dairong Li, Hong Wang, Yujuan He, Ju Cao, Yibing Yin
Infect Immun. 2011 February; 79(2): 867–878. Published online 2010 November 22.
doi: 10.1128/IAI.00717-10
PMCID: PMC3028838

161: c-Abl Phosphorylation of Mdm2 Facilitates Mdm2-Mdmx Complex Formation
David L. Waning, Jason A. Lehman, Christopher N. Batuello, Lindsey D. Mayo
J Biol Chem. 2011 January 7; 286(1): 216–222. Published online 2010 November 16.
doi: 10.1074/jbc.M110.183012

PMCID: PMC3012977

162: The scaffolding protein EBP50 regulates microvillar assembly in a phosphorylation-dependent manner

Damien Garbett, David P. LaLonde, Anthony Bretscher

J Cell Biol. 2010 October 18; 191(2): 397–413. doi: 10.1083/jcb.201004115

PMCID: PMC2958488

163: Crystal Structure of Cardiac-specific Histone Methyltransferase SmyD1 Reveals Unusual Active Site Architecture

Nualpun Sirinupong, Joseph Brunzelle, Jun Ye, Ali Pirzada, Lindsey Nico, Zhe Yang

J Biol Chem. 2010 December 24; 285(52): 40635–40644. Published online 2010 October 12.

doi: 10.1074/jbc.M110.168187

PMCID: PMC3003362

164: Isolation of SPINK6 in Human Skin: SELECTIVE INHIBITOR OF KALLIKREIN-RELATED PEPTIDASES

Ulf Meyer-Hoffert, Zhihong Wu, Tomasz Kantyka, Jan Fischer, Ties Latendorf, Britta

Hansmann, Joachim Bartels, Yinghong He, Regine Gläser, Jens-Michael Schröder

J Biol Chem. 2010 October 15; 285(42): 32174–32181. Published online 2010 July 28.

doi: 10.1074/jbc.M109.091850

PMCID: PMC2952218

165: The structure of FANCL, the catalytic subunit of the Fanconi Anemia core complex
Ambrose R. Cole, Laurence P.C. Lewis, Helen Walden

Nat Struct Mol Biol. Author manuscript; available in PMC 2010 September 1. Published in final

edited form as: Nat Struct Mol Biol. 2010 March; 17(3): 294–298. Published online 2010

February 14. doi: 10.1038/nsmb.1759

PMCID: PMC2929457

166: A Gate-Latch-Lock Mechanism for Hormone Signaling by Abscisic Acid Receptors

Karsten Melcher, Ley-Moy Ng, X. Edward Zhou, Fen-Fen Soon, Yong Xu, Kelly M. Suino-

Powell, Sang-Youl Park, Joshua J. Weiner, Hiroaki Fujii, Viswanathan Chinnusamy, Amanda

Kovach, Jun Li, Yonghong Wang, Jiayang Li, Francis C. Peterson, Davin R. Jensen, Eu-Leong

Yong, Brian F. Volkman, Sean R. Cutler, Jian-Kang Zhu, H. Eric Xu

Nature. Author manuscript; available in PMC 2010 June 3. Published in final edited form as:

Nature. 2009 December 3; 462(7273): 602–608. doi: 10.1038/nature08613

PMCID: PMC2810868

167: Pivotal Role of the C-terminal DW-motif in Mediating Inhibition of Pyruvate
Dehydrogenase Kinase 2 by Dichloroacetate

Jun Li, Masato Kato, David T. Chuang

J Biol Chem. 2009 December 4; 284(49): 34458–34467. Published online 2009 October 15.

doi: 10.1074/jbc.M109.065557

PMCID: PMC2797213

168: Distinct Amino Termini of Two Human HCS Isoforms Influence Biotin Acceptor Substrate Recognition

Maria Ingaramo, Dorothy Beckett

J Biol Chem. 2009 November 6; 284(45): 30862–30870. Published online 2009 September 9.

doi: 10.1074/jbc.M109.046201

PMCID: PMC2781485

169: Structural Studies of Soybean Calmodulin Isoform 4 Bound to the Calmodulin-binding Domain of Tobacco Mitogen-activated Protein Kinase Phosphatase-1 Provide Insights into a Sequential Target Binding Mode

Hiroaki Ishida, Mario Rainaldi, Hans J. Vogel

J Biol Chem. 2009 October 9; 284(41): 28292–28305. Published online 2009 August 10.

doi: 10.1074/jbc.M109.025080

PMCID: PMC2788880

170: Stabilization of Peptide Vesicles by Introducing Inter-Peptide Disulfide Bonds

Albert J. van Hell, Daan J. A. Crommelin, Wim E. Hennink, Enrico Mastrobattista

Pharm Res. 2009 September; 26(9): 2186–2193. Published online 2009 July 7.

doi: 10.1007/s11095-009-9933-z

PMCID: PMC2719749

171: Cdc7p-Dbf4p Regulates Mitotic Exit by Inhibiting Polo Kinase

Charles T. Miller, Carrie Gabrielse, Ying-Chou Chen, Michael Weinreich

PLoS Genet. 2009 May; 5(5): e1000498. Published online 2009 May 29.

doi: 10.1371/journal.pgen.1000498

PMCID: PMC2682205

172: Characterization of purified Sindbis Virus nsP4 RNA-dependent RNA Polymerase activity in vitro

Jon K. Rubach, Brian R. Wasik, Jonathan C. Rupp, Richard J. Kuhn, Richard W. Hardy, Janet L. Smith

Virology. Author manuscript; available in PMC 2011 June 3. Published in final edited form as:

Virology. 2009 February 5; 384(1): 201–208. Published online 2008 November 25.

doi: 10.1016/j.virol.2008.10.030

PMCID: PMC3107704

173: A DNA polymerase epsilon mutant that specifically causes +1 frameshift mutations within homonucleotide runs in yeast.

J M Kirchner, H Tran, M A Resnick

Genetics. 2000 August; 155(4): 1623–1632.

PMCID: PMC1461198

174: Norovirus Proteinase-Polymerase and Polymerase Are Both Active Forms of RNA-Dependent RNA Polymerase

Gaël Belliot, Stanislav V. Sosnovtsev, Kyeong-Ok Chang, Vijay Babu, Uzo Uche, Jamie J.

Arnold, Craig E. Cameron, Kim Y. Green

J Virol. 2005 February; 79(4): 2393–2403. doi: 10.1128/JVI.79.4.2393-2403.2005
PMCID: PMC546540

175: ClpS modulates but is not essential for bacterial N-end rule degradation
Kevin H. Wang, Robert T. Sauer, Tania A. Baker
Genes Dev. 2007 February 15; 21(4): 403–408. doi: 10.1101/gad.1511907
PMCID: PMC1804329

176: Genetic factors affecting the impact of DNA polymerase delta proofreading activity on
mutation avoidance in yeast.
H T Tran, N P Degtyareva, D A Gordenin, M A Resnick
Genetics. 1999 May; 152(1): 47–59.
PMCID: PMC1460598

177: Ni W, Xu SL, Tepperman JM, Stanley DJ, Maltby DA, Gross JD, Burlingame AL, Wang
ZY, Quail PH. A mutually assured destruction mechanism attenuates light signaling in
Arabidopsis. Science. 2014 Jun 6;344(6188):1160-4. doi:10.1126/science.1250778. PubMed
PMID: 24904166; PubMed Central PMCID: PMC4414656.

178: Fradet-Turcotte A, Canny MD, Escribano-Díaz C, Orthwein A, Leung CC, Huang H,
Landry MC, Kitevski-LeBlanc J, Noordermeer SM, Sicheri F, Durocher D. 53BP1 is a reader of
the DNA-damage-induced H2A Lys 15 ubiquitin mark. Nature. 2013 Jul 4;499(7456):50-4. doi:
10.1038/nature12318. Epub 2013 Jun 12. PubMed PMID:23760478; PubMed Central PMCID:
PMC3955401.

179: Bozza WP, Liang Q, Gong P, Zhuang Z. Transient kinetic analysis of USP2-catalyzed
deubiquitination reveals a conformational rearrangement in the K48-linked diubiquitin substrate.
Biochemistry. 2012 Dec 18;51(50):10075-86. doi:10.1021/bi3009104. Epub 2012 Dec 4.
PubMed PMID: 23211065; PubMed Central PMCID:

180: Hjerpe, R., Aillet, F., Lopitz-Otsoa, F., Lang, V., England, P., and Rodrigues, M.S. (2009)
Efficient protection and isolation of ubiquitylated proteins using tandem ubiquitin-binding
entities. EMBO Rep. 10,1250-1258.