



REFERENCES from peer reviewed publications

Imaging

Schröder, J., Benink, H., Dyba, M. and Los, G.V. (2009) In vivo labeling method using a genetic construct for nanoscale resolution microscopy. *Biophys J.* **96**, L1-3.

Johnsson, K. (2009) Visualizing biochemical activities in living cells. *Nat. Chem. Biol.*, **5**, 63–5.

Porter, F.W. and Palmenberg, A.C. (2009) Leader-induced phosphorylation of nucleoporins correlates with nuclear trafficking inhibition by cardioviruses. *J. Virol.* **83**, 1941–51.

Yamaguchi, K. *et al.* (2009) Pulse-chase experiment for the analysis of protein stability in cultured mammalian cells by covalent fluorescent labeling of fusion proteins. *Methods in Molecular Biology*, (in press).

Svendsen, S. *et al.* (2008) Spatial separation and bidirectional trafficking of proteins using a multi-functional reporter. *BMC Cell Biology*, **9**, 17 (<http://www.biomedcentral.com/1471-2121/9/17>).

Cao, Z. *et al.* (2008) Use of fluorescence-activated vesicle sorting for isolation of naked2-associated, basolaterally-targeted exocytic vesicles for proteomic analysis. *Mol. Cell Proteomics* **7**, 1651–67.

Yamaguchi, K., Ohara, O., Ando, A. and Nagase, T. (2008) Smurf1 directly targets hPEM-2, a GEF for Cdc42, via a novel combination of protein interaction modules in the ubiquitin-proteasome pathway. *Biological Chemistry* **389**, 405–13.

Cho, C., Reck-Peterson, S.L. and Vale, R.D. (2008) Regulatory ATPase sites of cytoplasmic dynein affect processivity and force generation. *J. Biol. Chem.* **283**, 25839–45.

Lin, M.Z. and Wang, L. (2008) Selective labeling of proteins with chemical probes in living cells. *Physiology* **23**, 131–41.

So, M.K., Yao, H. and Rao, J. (2008) HaloTag Protein-mediated specific labeling of living cells with quantum dots. *Biochemical and Biophysical Research Communications* **374**, 419.

Tanematsu, K. *et al.* (2008) Identification of ubiquitin ligase activity of RBCK1 and its inhibition by splice variant RBCK2 and protein kinase C β . *J. Biol. Chem.* **283**, 11575–85.

Wang, L. *et al.* (2008) The chaperone activity of heat shock protein 90 is critical for maintaining the stability of leucine-rich repeat kinase 2. *J. Neurosci.* **28**, 3384–91.

Hirakawa, Y. *et al.* (2008) Non-cell autonomous control of vascular stem cell fate by a CLE peptide/receptor system. *PNAS* **105**, 15208–13.

Beltrami, A. *et al.* (2008) Citrullination-dependent differential presentation of a self-peptide by HLA-B27 subtypes. *J. Biol. Chem.* **283**, 27189–99.



REFERENCES from peer reviewed publications

Ogawa, M., Shinohara, H., Sakagami, Y. and Matsubayashi, Y. (2008) Arabidopsis CLV3 peptide directly binds CLV1 ectodomain. *Science* **319**, 274.

Gautier, A. *et al.* (2008) An engineered protein tag for multiprotein labeling in living cells. *Chem. Biol.* **15**, 128–36.

Straszewski-Chavez, S.L. *et al.* (2007) XAF1 mediates tumor necrosis factor- α -induced apoptosis and X-linked inhibitor of apoptosis cleavage by acting through the mitochondrial pathway. *J. Biol. Chem.* **282**, 13059–72.

Esposito, A., Dohm, C. P., Bahr, M. and Wouters, F.S. (2007) Unsupervised fluorescence lifetime imaging microscopy for high content and high throughput screening. *Mol. Cell Proteomics* **6**, 1446–54.

Gerlach, R.G., Holzer, S.U., Jackel, D. and Hensel, M. (2007) Rapid engineering of bacterial reporter gene fusions by using Red recombination. *Appl. Environ. Microbiol.* **73**, 4234–42.

Honma, N. (2007) HaloTag: a novel technology for in vivo bioimaging and protein functional analysis. *Tanpakushitsu Kakusan Koso.* **52**, 1575–80.

Zhang, Y. *et al.* (2006) HaloTag protein-mediated site-specific conjugation of bioluminescent proteins to quantum dots. *Angewandte Chemie International Edition.* **45**, 4936–40.

Lang, C., Schulze, J., Mendel, R.R. and Hansch, R. (2006) HaloTag®: a new versatile reporter gene system in plant cells. *Journal of Experimental Botany.* **B**, 2985–92.

Reck-Peterson, S.L. *et al.* (2006) Single-molecule analysis of dynein processivity and stepping behavior. *Cell* **126**, 335–48.

Protein Immobilization

Nath, N. *et al.* (2008) Improving protein array performance: focus on washing and storage conditions. *Journal of Proteome Research* **7 (10)**, 4475–82.

Hata, T. and Nakayam, M. (2007) Rapid single-tube method for small-scale affinity purification of polyclonal antibodies using HaloTag® Technology. *J. Biochemical and Biophysical Methods* **70**, 679–82.

Shinohara, H. and Matsubayashi, Y. (2007) Functional immobilization of plant receptor-like kinase onto microbeads towards receptor array construction and receptor-based ligand fishing. *The Plant Journal* **52(1)**:175–84.

Additional Citations

Kardon, J.R., Reck-Peterson, S.L. and Vale, R.D. (2009) Regulation of the processivity and intracellular localization of *Saccharomyces cerevisiae* dynein by dynactin. *PNAS* **106**, 5669–74.

Los, G.V. *et al.* (2008) HaloTag®: a novel protein labeling technology for cell imaging and protein analysis. *ACS Chemical Biology* **3**, 373–82.



REFERENCES from peer reviewed publications

Nagase, T. *et al.* (2008) Exploration of human ORFeome: high-throughput preparation of ORF clones and efficient characterization of their protein products. *DNA Research* **15**, 137–49.

Ogawa, M., Shinohara, H., Sakagami, Y. and Matsubayashi, Y. (2008) Arabidopsis CLV3 peptide directly binds CLV1 ectodomain. *Science* **319**, 294.

Lee, J., Yu, P., Xiao, X. and Kodadek, T. (2008) A general system for evaluating the efficiency of chromophore-assisted light inactivation (CALI) of proteins reveals Ru(II) tris-bipyridyl as an unusually efficient warhead. *Molecular Biosys.* **4**, 59–65.

Los, G.V. and Wood, K. (2007) The HaloTag®: a novel technology for cell imaging and protein analysis. *Methods Mol Biol.* **356**, 195–208.

Temple, G. *et al.* (2006) From genome to proteome: developing expression clone resources for the human genome. *Hum. Mol. Genet.* **15**, R31–43.

