

Publications using the in situ Proximity Ligation Assay Technology in order of  
publication date

Spears M, Bartlett JM. Human Epidermal Growth Factor Receptor Dimerization Analysis in Breast Cancer Diagnosis: Potential for Improving Testing Accuracy and Treatment Selection. *Mol Diagn Ther*. 2009 Dec 1;13(6):359-365.PMID: 19925033

Kii I, Incorporation of tenascin-C into the extracellular matrix by periostin underlies an extracellular meshwork architecture. *J Biol Chem*. 2009 Nov 3.

Hasumi Y, Baba M, Ajima R, Hasumi H, Valera VA, Klein ME, Haines DC, Merino MJ, Hong SB, Yamaguchi TP, Schmidt LS, Linehan WM. Homozygous loss of BHD causes early embryonic lethality and kidney tumor development with activation of mTORC1 and mTORC2. *Proc Natl Acad Sci U S A*. 2009 Nov 3;106(44):18722-7.

Leuchowius KJ, Jarvius M, Wickström M, Rickardson L, Landegren U, Larsson R, Söderberg O, Fryknäs M, Jarvius J. High content screening for inhibitors of protein interactions and post-translational modifications in primary cells by proximity ligation. *Mol Cell Proteomics*. 2009 Oct 27.

Hervouet E, Dnmt3/transcription factor interactions as crucial players in targeted DNA methylation. *Epigenetics*. 2009 Oct 21;4(7).

Robinson SD, Alphavbeta3-integrin limits the contribution of neuropilin-1 to VEGF-induced angiogenesis. *J Biol Chem*. 2009 Oct 16.

Smith R, Mutant huntingtin interacts with {beta}-tubulin and disrupts vesicular transport and insulin secretion. *Hum Mol Genet*. 2009 Oct 15;18(20):3942-54.

Leuchowius KJ, Weibrecht I, Landegren U, Gedda L, Söderberg O. Flow cytometric in situ proximity ligation analyses of protein interactions and post-translational modification of the epidermal growth factor receptor family. *Cytometry A*. 2009 Oct;75(10):833-9

Tim Conze, Ana Sofia Carvalho, Ulf Landegren, Raquel Almeida, Celso A. Reis, Leonor David, and Ola Söderberg  
MUC2 mucin is a major carrier of the cancer-associated sialyl-Tn antigen in intestinal metaplasia and gastric carcinomas.. *Glycobiology*, Advance Access published on October 8, 2009

Per Svenningsen, Torben R. Uhrenholt, Yaseelan Palarasah, Karsten Skjoedt, Boye L. Jensen, and Ole Skott.  
Prostasin-dependent activation of epithelial Na<sup>+</sup> channels by low plasmin concentrations. *Am J Physiol Regul Integr Comp Physiol* 2009 Sept 30

Koos B, Paulsson J, Jarvius M, Sanchez BC, Wrede B, Mertsch S, Jeibmann A, Kruse A, Peters O, Wolff JE, Galla HJ, Söderberg O, Paulus W, Ostman A, Hasselblatt M.

Platelet-Derived Growth Factor Receptor Expression and Activation in Choroid Plexus Tumors. **Am J Pathol**. 2009 Aug 28. [Epub ahead of print]

Yadavilli S, Mayo LD, Higgins M, Lain S, Hegde V, Deutsch WA.  
Ribosomal protein S3: A multi-functional protein that interacts with both p53 and MDM2 through its KH domain.**DNA Repair** (Amst). 2009 Aug 3. Epub ahead of print

Karl-Johan Leuchowius, Irene Weibrecht, Ulf Landegren, Lars Gedda, Ola Söderberg.  
Flow cytometric in situ proximity ligation analyses of protein interactions and post-translational modification of the epidermal growth factor receptor family. **Cytometry Part A**. Published Online 31 Jul 2009.

Yadav AK, Renfrow JJ, Scholtens DM, Xie H, Duran GE, Bredel C, Vogel H, Chandler JP, Chakravarti A, Robe PA, Das S, Scheck AC, Kessler JA, Soares MB, Sikic BI, Harsh GR, Bredel M.

Monosomy of chromosome 10 associated with dysregulation of epidermal growth factor signaling in glioblastomas. **JAMA**. 2009 Jul 15;302(3): 276-89.

Tuomi S, Mai A, Nevo J, Laine JO, Vilkki V, Ohman TJ, Gahmberg CG, Parker PJ, Ivaska

PKCepsilon regulation of an alpha5 integrin-ZO-1 complex controls lamellae formation in migrating cancer cells.**J.Sci Signal**. 2009 Jun 30;2(77):ra32.PMID: 19567915 [PubMed - in process]

Massinen S, Tammimies K, Tapia-Páez I, Matsson H, Hokkanen ME, Söderberg O, Landegren U, Castrén E, Gustafsson JA, Treuter E, Kere J.

Functional interaction of DYX1C1 with estrogen receptors suggests involvement of hormonal pathways in dyslexia. **Hum Mol Genet**. 2009 May 7. [Epub ahead of print]

Tan SH, Pal M, Tan MJ, Wong MH, Tam FU, Teo JW, Chong HC, Tan CK, Goh YY, Tang MB, Cheung PC, Tan NS.

Regulation of cell proliferation and migration by TAK1 via transcriptional control of von hippel-lindau tumor suppressor. **J Biol Chem**. 2009 May 6. [Epub ahead of print]

E Laane, K Pokrovskaja Tamm, E Buentke, K Ito, P Khahariza, J Oscarsson, M Corcoran, A-C Björklund, K Hultenby, J Lundin, M Heyman, S Söderhäll, J Mazur, A Porwit, PP Pandolfi, B Zhivotovsky, T Panaretakis, and D Grandér.

Cell death induced by dexamethasone in lymphoid leukemia is mediated through initiation of autophagy. Cell Death and Differentiation advance online publication, 24 April 2009; doi10.1038/cdd.2009.46

Yamazaki T, Yoshimatsu Y, Morishita Y, Miyazono K, Watabe T.

COUP-TFII regulates the functions of Prox1 in lymphatic endothelial cells through direct interaction. *Genes Cells*. 2009 Mar;14(3):425-34. Epub 2009 Feb 4.

Mellberg S, Dimberg A, Bahram F, Hayashi M, Rennel E, Ameer A, Westholm JO, Larsson E, Lindahl P, Cross MJ, Claesson-Welsh L.

Transcriptional profiling reveals a critical role for tyrosine phosphatase VE-PTP in regulation of VEGFR2 activity and endothelial cell morphogenesis. *FASEB J.* 2009 Jan 9. [Epub ahead of print] PMID: 19136612

Greenberg JI, Shields DJ, Barillas SG, Acevedo LM, Murphy E, Huang J, Scheppke L, Stockmann C, Johnson RS, Angle N, Cheresh DA.

A role for VEGF as a negative regulator of pericyte function and vessel maturation. *Nature.* 2008 Nov 9. [Epub ahead of print]

Taro Matsumoto, Petter Schiller, Lothar C. Dieterich, Fuad Bahram, Yuji Iribe, Ulf Hellman, Charlotte Wikner, Gordon Chan, Lena Claesson-Welsh and Anna Dimberg. Ninein is expressed in the cytoplasm of angiogenic tip-cells and regulates tubular morphogenesis of endothelial cells. *Arteriosclerosis, Thrombosis, and Vascular Biology.* Published online Sep 4, 2008

Vermeer MH, van Doorn R, Dijkman R, Mao X, Whittaker S, van Voorst Vader PC, Gerritsen MJ, Geerts ML, Gellrich S, Söderberg O, Leuchowius KJ, Landegren U, Out-Luiting JJ, Knijnenburg J, Ijszenga M, Szuhai K, Willemze R, Tensen CP.

Novel and highly recurrent chromosomal alterations in Sézary syndrome. *Cancer Res.* 2008 Apr 15

Szabo T., Vanderheyden V., Parys JB., De Smedt H., Rietdorf K., Kotelevets L., Chastre E., Khan F., Landegren U., Söderberg O., Bootman MD., Roderick HL.

Phosphorylation of inositol 1,4,5-trisphosphate receptors by protein kinase B/Akt inhibits Ca<sup>2+</sup> release and apoptosis. *Proc Natl Acad Sci U S A.* 2008 Feb 19;105(7):2427-32. Epub 2008 Feb 4.

Jarvis, M., Paulsson, J., Weibrecht, I., Leuchowius, K., Andersson, A-C., Wahlby, C., Gullberg, M., Sjöblom, T., Markova, B., Ostman, A., Landegren, U., & Soderberg, O., In situ detection of phosphorylated PDGF receptor beta using a generalized proximity ligation method. *Mol. Cell Proteomics*, 2007

Soderberg, O., Gullberg, M., Jarvis, M., Ridderstrale, K., Leuchowius, K. J., Jarvis, J., Wester, K., Hydbring, P., Bahram, F., Larsson, L. G. & Landegren, U.

Direct observation of individual endogenous protein complexes in situ by proximity ligation. *Nat. Methods*, 2006.

Jarvis, J., Melin, J., Goransson, J., Stenberg, J., Fredriksson, S., Gonzalez-Rey, C., Bertilsson, S. & Nilsson, M.,

Digital quantification using amplified single-molecule detection. *Nat. Methods*, 2006, 3, 725-7.