

## January 2012 Webinar: Overcoming Challenges of Protein Analysis in Mammalian Systems

### Key Websites Cited:

#### Kazusa DNA Research Institute

Human ORFs N-terminal fusion constructs in HaloTag® Flexi® vector pFN21A for expression in mammalian cells

<http://www.kazusa.or.jp/kop/dsearch-e//>

#### GeneCopoeia

Human and mouse ORFs in N- /C- terminal fusion constructs available in OmicsLink™ vectors for expression in mammalian cells <http://www.genecopoeia.com/product/halo/>

### Publications Cited:

#### HaloTag-based purification of functional human kinases from mammalian cells.

RF Ohana, R Hurst, J Vidugiriene, MR Slater, KV Wood, and M Urh  
(2011) *Protein Expr Purif*, **76**(2): 154-64.

<http://www.sciencedirect.com/science/article/pii/S1046592810003244>

#### Spatial separation and bidirectional trafficking of proteins using a multi-functional reporter.

S Svendsen, C Zimprich, MG McDougall, DH Klaubert, and GV Los  
(2008) *BMC Cell Biol*, **9**: 17.

<http://www.biomedcentral.com/1471-2121/9/17>

#### A functional analysis of the CREB signaling pathway using HaloCHIP-chip and high throughput reporter assays.

DD Hartzell, ND Trinklein, J Mendez, N Murphy, SF Aldred, K Wood, and M Urh  
(2009) *BMC Genomics*, **10**: 497.

<http://www.biomedcentral.com/1471-2164/10/497>

#### Examining the Complexity of Human RNA Polymerase Complexes using HaloTag Technology Coupled to Label Free Quantitative Proteomics.

Daniels, D.L., *et al.*

(2012) *J Proteome Res*, Article ASAP

<http://pubs.acs.org/doi/full/10.1021/pr200459c?prevSearch=daniels%252C%2BD&searchHistoryKey=>

#### Structure of the 70S Ribosome Complexed with mRNA and tRNA.

Maria Selmer, Christine M. Dunham, Frank V. Murphy, IV, Albert Weixlbaumer, Sabine Petry, Ann C. Kelley, John R. Weir, and V. Ramakrishnan

(2006) *Science*, **313**: 1935-1942.

<http://www.sciencemag.org/content/313/5795/1935.full.pdf>