



V3 WESTERN WORKFLOW™

## Housekeeping Protein (HKP) Publications

### Research Papers on the Pitfalls of Using HKP as Loading Controls in Western Blots

**Avoiding pitfalls of internal controls: validation of reference genes for analysis by qRT-PCR and western blot throughout rat retinal development.**

Rocha-Martins M, Njaine B, Silveira MS (2012). PLoS One 7, e43028.

**Using soybean trypsin inhibitor as an external loading control for western blot analysis of tear proteins: application to corneal disease.**

You J, Hodge C, Wen L, McAvoy JW, Madigan MC, Sutton G (2012). Exp Eye Res 99, 55–62.

**Stability of reference proteins in human placenta: general protein stains are the benchmark.**

Lanoix D, St-Pierre J, Lacasse AA, Viau M, Lafond J, Vaillancourt C (2012). Placenta 33, 151–156.

**Uncovering suitable reference proteins for expression studies in human adipose tissue with relevance to obesity.**

Pérez-Pérez R, López JA, García-Santos E, Camafeita E, Gómez-Serrano M, Ortega-Delgado FJ, Ricart W, Fernández-Real JM, Peral B (2012). PLoS One 7, e30326.

**Identification and validation of rice reference proteins for western blotting.**

Li X, Bai H, Wang X, Li L, Cao Y, Wei J, Liu Y, Liu L, Gong X, Wu L, Liu S, Liu G (2011). J Exp Bot 62, 4763–4772.

**Housekeeping genes: expression levels may change with density of cultured cells.**

Greer S, Honeywell R, Geletu M, Arulanandam R, Raptis L (2010). J Immunol Methods 355, 76–79.

**Absence of GAPDH regulation in tumor cells of different origin under hypoxic conditions in vitro.**

Said HM, Polat B, Hagemann C, Anacker J, Flentje M, Vordermark D (2009). BMC Res Notes 2, 8.

**Expression of four housekeeping proteins in elderly patients with schizophrenia.**

Bauer DE, Haroutunian V, McCullumsmith RE, Meador-Woodruff JH (2009). J Neural Transm 116, 487–491.

**Beta-tubulin is a more suitable internal control than beta-actin in western blot analysis of spinal cord tissues after traumatic injury.**

Liu NK and Xu XM (2006). J Neurotrauma 12, 1794–1801.

**Beta-actin is not a reliable loading control in western blot analysis.**

Dittmer A and Dittmer J (2006). Electrophoresis 27, 2844–2845.

**Housekeeping proteins: a preliminary study illustrating some limitations as useful references in protein expression studies.**

Ferguson RE, Carroll HP, Harris A, Maher ER, Selby PJ, Banks RE (2005). Proteomics 5, 566–571.

**Glyceraldehyde-3-phosphate dehydrogenase varies with age in glycolytic muscles of rats.**

Lowe DA, Degens H, Chen KD, Alway SE (2000). J Gerontol A Biol Sci Med Sci 55, B160–164.

**BIO-RAD**