

TnT[®] *in vitro* Transcription/ Translation Systems

One system, multiple applications with TnT[®] *in vitro* Transcription/Translation Systems

Promega TnT[®] Coupled Transcription/Translation Systems offer a one step, easy to perform protocol with significant time savings over cellular approaches to protein expression. Proteins expressed using these systems can be used for a variety of applications relating to their relationship with other molecules and function.

References

Protein Pull-downs

- Ma, W. *et al.* (2010) *J. Virol.* **84**, 2122-33.
- Kim, S.M. *et al.* (2010) *Nucl. Acid. Res.* **38**, 6389-403.
- Camacho, L. *et al.* (2009) *J. Cell. Sci.* **122**, 4383-92.
- Houshmandi, S. *et al.* (2009) *Mol. Cell Biol.* **29**, 1472-86.

Gel Shift (EMSA)

- Buas, M. *et al.* (2010) *J. Biol. Chem.* **285**, 1249-58.
- Lin, Y. *et al.* (2009) *Drug. Metab. Dispos.* **37**, 1295-304.
- Bollig, F. *et al.* (2009) *Develop.* **136**, 2883-92.

Chloroplast Import Assays

- Tanz, S. *et al.* (2009) *Plant. Phy.* **150**, 1515-29.

Co-immunoprecipitation

- Xi, Gang. *et al.* (2010) *J. Biol. Chem.* **285**, 6937-51.
- Martin, L. and Tremblay, J. (2009) *J. Mol. Endocrinol.* **42**, 119-29.
- Makarenkova, H. *et al.* (2009) *J. Biol. Chem.* **284**, 14866-74.
- Marchetti, B. *et al.* (2009) *J. Gen. Virol.* **90**, 2865-70.

Ubiquitination Assays

- Cen, B. *et al.* (2010) *J. Biol. Chem.* **285**, 29128-37.
- Zhong, B. *et al.* (2010) *J. Immunol.* **184**, 6249-55.
- Fu, W. *et al.* (2009) *J. Biol. Chem.* **284**, 13987-14000.

Protein Folding

- Berry, J-L and Bulleid, N. (2009) *BioSci. Horz.* **2**, 13-21.

Sumoylation Assays

- Rytinki, M. *et al.* (2009) *J. Biol. Chem.* **284**, 26184-93.
- Klein, U. *et al.* (2009) *Mol. Cell. Biol.* **20**, 410-18.
- Seo, W. and Ziltener, H. (2009) *Blood*, **17**, 3567-77.

Inhibition of Translation

- Zhang, Y. and Inouye, M. (2009) *J. Biol. Chem.* **284**, 6627-38.
- Liu, Y. *et al.* (2009) *J. Biol. Chem.* **284**, 5859-68.

Membrane Association

- Bintintan, I. and Meyers, G. (2010) *J. Biol. Chem.* **285**, 8572-84.
- Al, L-S. *et al.* (2009) *J. Virol.* **83**, 9923-39.
- Stapleford, K. *et al.* (2009) *J. Virol.* **83**, 4498-507.