



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

***PTP PEST*** (Human, recombinant, N-terminal GST tag)  
Catalog #: 30022

**Description:** Protein tyrosine phosphatase, Catalytic domain, (a.a. 2-300) (Gene bank accession number NM\_002835) with N-terminal GST tag, MW=61.4 kDa, expressed in an E.coli expression system.

**Source:** E.Coli

**Formulated in** 25 mM Tris-HCl, pH 8.0, 75 mM NaCl, 0.05% Tween-20, 50% glycerol, 2 mM EDTA, 1 mM DTT, 10 mM glutathione.

**Stability:** >6 months at -80°C

**Purity:** > 90% by SDS-PAGE

**Specific Activity:** 1 U/ug. One unit will hydrolyze 1 nmol p-nitrophenyl phosphate per minute at pH 7.4 and 30°C. Assay buffer: 50 mM HEPES, pH 7.4, 2 mM EDTA, 3 mM DTT, 100 mM NaCl, 50 mM pNPP.

**Application:** Useful for the study of enzyme kinetics, regulation, to dephosphorylate target substrates and for screening inhibitors.

### **References:**

1. J.F. Cote, et al., *J Biol Chem.* 2002, 277: 2973-86.
2. K. Horsch et al., *Mol Endocrinol.* 2001, 12: 2182-96.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)