



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

***PTP1B*** (rat, recombinant, N-terminal GST tag)  
Catalog #: 30011

**Description:** Protein tyrosine phosphatase, catalytic domain (a.a. 1-321), with N-terminal GST tag, MW=63.3 kDa, expressed in an E.coli expression system.

**Source:** E.Coli

**Formulation:** Recombinant proteins were eluted in 25 mM Tris-HCL, 75 mM NaCl, pH 8.0, 0.05% Tween, 10 mM glutathione, 1 mM DTT and 50% glycerol. Liquid at – 20°C.

**Stability:** >6 months at -80°C

**Purity:** > 95% by SDS-PAGE

**Specific Activity:** The specific activity of rat PTP1B was determined using pNPP.  
Enzyme reaction condition: 20 mM pNPP, 5 min incubation at 30°C, 2 ug/ml enzyme.

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

### **References:**

1. Frangioni et al., *Cell*, 1992, 68: 545
2. M. Elchebly et al., *Science* 1999, 283: 1544

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)