



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### ***PTP1B (mouse)***

**Catalog #:** 30012

**Lot#:** 1000    **Conc.:** 1.5 mg/ml

**Expression System:** Murine recombinant human PTP1B Protein tyrosine phosphatase, also known as PTPN1, catalytic domain, (a.a. 1-321), (Genbank accession no. NM\_011201.3) with an N-terminal GST-tag, MW=63.3 kDa, expressed in *E. coli*.

**Purity:** The purity is >95%. Purified using glutathione-agarose beads. The purified GST fusion proteins were analyzed on Coomassie-stained SDS gel.

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

**Activity:** The activity of mouse PTP1B was determined using pNPP.

**Assay Conditions:** 50 mM HEPES, pH 7.4, 2 mM EDTA, 3 mM DTT, 100 mM NaCl, 20 mM pNPP, 5 min incubation at 30°C, using 2 µg/ml enzyme

**Unit Definition:** One unit will hydrolyze 1 nmol p-nitrophenyl phosphate per minute at pH 7.4 and 30°C.

**Stability:** Stable for >6 months at -80°C from date of shipment.

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

#### **References:**

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

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
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)