



6044 Cornerstone Court West, Suite 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**

### ***PTP-Bas***

Catalog #: 30024  
Lot #: 2000 Conc.: 1.5 mg/ml

**Description:** Human, recombinant PTP-Bas, Catalytic domain, (a.a. 2091-2490), (Gene bank accession number NM\_080685) with N-terminal GST tag, MW=71.8 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:** >80% by SDS-PAGE.

**Specific Activity:** 5 U/μg. 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, and 100 mM NaCl. The specific activity of Bas was determined using pNPP. Enzyme reaction condition: 50 mM HEPES, pH 7.4, 2 mM EDTA, 3 mM DTT, 100 mM NaCl, 20 mM pNPP, and 20 ng/μl enzyme, for 5 min at 30°C.

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

#### **References:**

1. W. Kimber, *et al.*, *Biochem. J.* 2003, **376**: 525-535
2. N. Kachel, *et al.* *J.Mol. Biol.* 2003, **334**: 143-155

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)