



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

Data Sheet

PTP1B (human, recombinant, N-terminal GST tag)
Catalog #: 30010

Description: Protein tyrosine phosphatase, catalytic domain, (a.a. 1-321), (Gene bank accession no. M31724) with N-terminal GST tag, MW=63.3 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: > 95% by SDS-PAGE

Specific Activity: 20 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. 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
Please visit our website at: www.bpsbioscience.com