



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 (mouse, recombinant, N-terminal GST tag)
Catalog #: 30012

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 mouse 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
Please visit our website at: www.bpsbioscience.com