



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

PRL-3 (human, recombinant, N-terminal GST tag)
Catalog #: 30002

Description: Protein tyrosine phosphatase, catalytic domain, phosphatase of regenerating liver3. Catalytic domain, (a.a. 2-173), (Gene bank accession number NM_032611) with N-terminal GST tag, MW=45.4 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: >3 months at -80°C

Purity: > 95% by SDS-PAGE

Specific Activity: 0.9 U/ug. One unit will hydrolyze 1 pmol 6, 8-difluoro-4-methyl umbelliferyl phosphate (DiFMUP) per minute at pH 6.3 and 30°C. Assay buffer: 50 mM Bis-Tris, pH 6.3, 2 mM EDTA, 2 mM DTT, 100 mM DiFMUP.

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

References:

1. S. Saha, et al., *Science* 2001, 294: 1343.
2. U.A. Miskad et al., *Pathobiology* 2004, 71: 176-84.

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