



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

DEP 1 (Human, recombinant, N-terminal GST tag)
Catalog #: 30050

Description: Protein tyrosine phosphatase. (density-enhanced phosphatase-1, PTP-eta, CD148), Catalytic domain, (a.a. 1024-1306) (Gene bank accession no. NM_U10886) with N-terminal GST tag, MW=58.9 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: 114 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, and 0.5 ug/ml enzyme.

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

References:

1. A. Massa, et al., *J Biol Chem.* 2004, 279: 29004-12
2. S. Kellie, et al., *J Cell Sci.* 2004, 117 :609-18

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