



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

Bas (Human, recombinant, N-terminal GST tag)
Catalog #: 30024

Description: Protein tyrosine phosphatase. Catalytic domain (aa 2091-2490) (Gene bank accession number NM_080685) with a GST-tag expressed in E.coli. Purified using glutathione-agarose beads.

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: > 80% by SDS-PAGE

Specific Activity: 5 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. The specific activity of Bas was determined using pNPP. Enzyme reaction condition: 20 mM pNPP, 5 min incubation at 30°C, 1 ug enzyme.

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

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