

Data Sheet

PTPβ

Human, recombinant, N-terminal GST tag
Catalog #: 30042
Lot #: 141117 **Conc.:** 4.8 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0.
 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 16
 mM glutathione and 20% glycerol

Stability: At least 6 months at -80°C .
 Avoid freeze/thaw cycles. Storing diluted
 enzyme is not recommended, if necessary,
 use carrier protein (BSA 0.1 – 0.5%).

References:

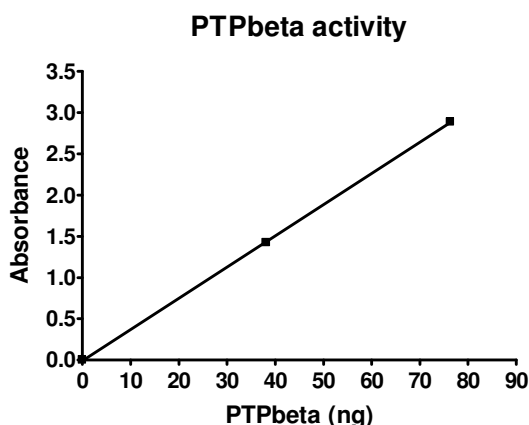
1. I. K. Lund et al., *J Biol Chem.* 2004,
279: 24226-35.
2. G.H. Peters et al., *Biochimie.* 2003,
85: 527-34.

Description: PTPβ (protein tyrosine
 phosphatase beta), catalytic domain, (a.a.
 1675-1996) with N-terminal GST tag,
 MW=63.7 kDa, expressed in an *E. coli*
 expression system.

Specific Activity: 55 nmole/min/μg.
 Assay conditions: 20 mM HEPES, pH 7.2,
 2 mM EDTA, 1 mM DTT, 25 mM pNPP
 and PTPβ. Reaction was done at room
 temperature for 10 minutes, and monitored
 at Abs 405 nm continuously. ϵ of pNPP =
 $18,000 \text{ M}^{-1}\text{cm}^{-1}$

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

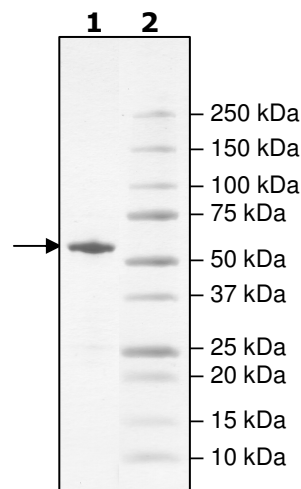
Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2.4 μg PTPbeta
Lane 2:
 Protein Marker

MW: 64 kDa
Purity: ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com