DATA SHEET





PTP 1B

Protein Tyrosine Phosphatase 1B human, recombinant, E. coli

Cat. No.	Amount
PR-931	20 μg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 37.3 kDa Accession number: NP_002818 Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM HEPES pH 7.4, 40 mM NaCl, 1 mM

EDTA, 1 mM DTT and 2 mM beta-Mercaptoethanol)

Description:

PTP 1B is the prototype nontransmembrane protein tyrosine phosphatase. It is a ubiquitous protein that was originally identified in human placenta. This highly active enzyme is well suited for the study of tyrosine phosphatase kinetics and substrate specificity, as well as for the screening of inhibitors.

1 nmol of PTP 1B will hydrolyze 37 nmol of pNPP per minute at 30°C.

Selected References:

Montalibet et al. (2004) Using yeast to screen for inhibitors of protein tyrosine phosphatase 1B. Biochem. Pharmacol. 68:1807.

Puius et al. (1997) Identification of a second aryl phosphatebinding site in protein-tyrosine phosphatase 1B: a paradigm for inhibitor design. Proc. Natl. Acad. Sci. USA. 94:13420.

Liu et al. (1996) Direct binding of the proline-rich region of protein tyrosine phosphatase 1B to the Src homology 3 domain of p130 (Cas). J. Biol. Chem. **271**:31290.

