

PTP 1B

Protein Tyrosine Phosphatase 1B

human, recombinant, *E. coli*

Cat. No.	Amount
PR-931	20 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50 mM HEPES pH 7.4, 40 mM NaCl, 1 mM EDTA and 1 mM DTT.

Activity

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

Purity

>95% by SDS-PAGE

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.

Selected References:

- 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.
- Puius *et al.* (1997) Identification of a second aryl phosphate binding site in protein-tyrosine phosphatase 1B: a paradigm for inhibitor design. *Proc. Natl. Acad. Sci. USA.* **94**:13420.
- Montalibet *et al.* (2004) Using yeast to screen for inhibitors of protein tyrosine phosphatase 1B. *Biochem. Pharmacol.* **68**:1807.