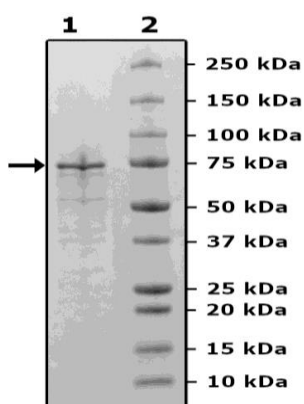


Product Information

Description	Protein tyrosine phosphatase nonreceptor, type 1 (PTP1B, also known as PTPN1), full length encompassing amino acids 1-435 (end). This construct contains an N-terminal GST-tag, linked to the protein via a short peptide corresponding to a thrombin cleavage site. The recombinant protein was affinity purified and tested for activity.
Species	Human
Construct:	PTP1B (GST-Th-1-435(end))
Concentration:	0.09 mg/ml
Expression System:	<i>E. coli</i>
Purity:	65%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
MW:	77 kDa
Genbank Accession:	M31724
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	2000 pmol/min/μg
Assay Conditions:	20 mM HEPES, pH 7.2, 2 mM EDTA, 1 mM DTT, 25 mM pNPP. The assay was done at room temperature for 1 hour and monitored at Abs ₄₁₅ nM. ϵ pNPP = 18,000 M ⁻¹ cm ⁻¹ with various amounts of PTP1B.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PTP1B Activity

