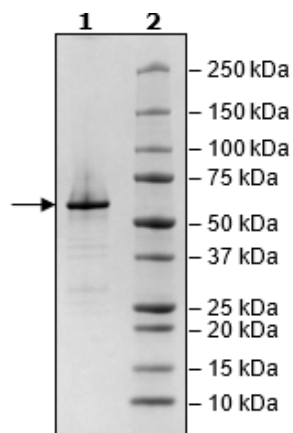


Product Information

Description:	Recombinant Mouse PTP1B (Protein tyrosine phosphatase), encompassing amino acids 1-321 (catalytic domain). This construct contains an N-terminal GST-tag. This recombinant protein was affinity purified.
Species	Mouse
Construct:	PTP1B (GST-1-321) (Mouse)
Concentration:	1.18 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	25 mM Tris-HCl, 75 mM NaCl, pH 8.0, 0.05% Tween-20, 10 mM glutathione, 1 mM DTT, and 50% glycerol
MW:	63 kDa
Genbank Accession:	NM_011201
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:	350 pmole/min/μg
Assay Conditions:	Assay was done according to PTP1B (Catalytic Domain) Colorimetric Assay Kit (BPS Bioscience #30019) 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



Mouse PTP1B Activity

