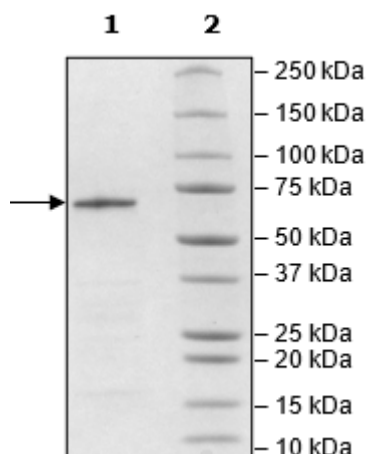


Product Information

Description:	Human recombinant SHP-2 (Src homology domain 2 containing tyrosine phosphatase, also known as protein tyrosine phosphatase N11, PTPN11), full length encompassing amino acids 2-593 (end). This construct contains an N-terminal His-tag (6xHis). The recombinant protein was affinity purified and tested for activity.
Species	Human
Construct:	SHP-2 (His-2-593(end))
Concentration:	0.18 mg/ml
Expression System:	<i>E. coli</i>
Purity:	83%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	70 kDa
Genbank Accession:	NM_002834
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:	≥50 pmol/min/μg
Assay Conditions:	Experiment was done according to SHP-2 (Full-Length) Homogeneous Assay Kit (BPS Bioscience #79330) with various amounts of SHP-2.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



SHP-2 Activity

