DATA SHEET





SHP-1^{GST} - full length

Src homology 2 domain Phosphatase-1 human, recombinant, *E. coli*

Cat. No.	Amount	
PR-947	20 µg	
	116.0 97.4 66.2	
	37.6	

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: 92 kDa

Accession number: NP_002822.2

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM Tris-HCl pH 8.0, 100 mM NaCl and 1 mM DTT)

Description:

The N-terminal GST-tagged fusion protein was expressed in E. coli and purified by affinity chromatography with GSH-beads. The GST-tag influences to some degree the stimulation by ligands of the N-terminal SH2-domain can dimerize and even be phosphorylated. The enzyme should only be used in diluted solutions or by adding 10% glycerol. SHP-1 (Src homology-2 containing protein tyrosine phosphatase-1) is a non-receptor protein tyrosine phosphatase with two phosphotyrosine binding domains. N- and C-terminal tandem SH2 domains lie N-terminal to the catalytic domain (PTP). In the unstimulated state interaction of the N-terminal SH2 domain with the catalytic domain leads to self inhibition. Natural ligand sequences from cytosolic parts of receptors, signal and scaffold proteins or synthetic phosphotyrosine peptides stimulate the phosphatase activity. Thus, SHP-1 acts as negative regulator in the signaling of various receptors, including erythropoietin receptor, IL3-receptor, CSF-1 receptor, B-cell receptor and c-Ros. SHP-1 prefers as substrate such proteins which are phosphorylated from the SRC-kinase. SHP-1 can act as tumor suppressor or can inhibit the processing of some immune cells.

Activity:

Basal activity (unstimulated): 130 pmol Pi/min/ μ g; Phosphatase activity was determined with pNPP as substrate at pH 7.5 and 30°C.(see PR-944) The enzyme was used in final concentrations of 15 and 30 nM.

Selected References:

Zhang *et al.* (2000) Roles of SHP-1 tyrosine phosphoatase in the negative regulation of Cell signallimg. *Immunology* **12**:361.

Keilhack *et al.* (2001) Negative regulation of Ros receptor tyrosine kinase signaling. An epithelial function of the SH2 domain protein tyrosine phosphatase SHP-1. *J. Cell Biol.* **152**:325.

Li et al. (2000) Form, function, and regulation of protein tyrosine phosphatases and their involvement in human diseases. Semin. Immunol. **12**:75.

Wu *et al.* (2003) SHP-1 suppresses cancer cell growth by promoting degradation of JAK kinases. *J. Cell Biochem.* **90**:1026.

Kamata *et al.* (2003) Src-homology 2 domain containing tyrosine phosphatase SHP-1 controls the development of allergic airway inflammation. *J. Clin. Invest.* **111**:109.

Yang et al. (1998) Crystal structure of the catalytic domain of protein-tyrosine phosphatase SHP-1. J. Biol. Chem. **273**:28199.

Yang et al.<7i> (2000) Structural basis for substrate specificity of protein-tyrosine phosphatase SHP-1. <i<J. Biol. Chem. **275**:4066.

Frank et al. (2004) Effective Dephosphorylation of Src Substrates by SHP-1. J. Biol. Chem. **279**:11375.



Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com