

# Shp-1 (human), (recombinant)

Shp-1 is one of two human "Src Homology-2 (SH2) containing phosphatases", enzymes in which two regulatory phosphotyrosine binding domains (N-SH2 & C-SH2) lie N-terminal to the catalytic (PTP) domain. Recruitment of Shp-1 (Src Homology-2 (SH2) containing phosphatase-1) to cell-surface receptors is mediated by SH2-domain binding to "Immuno-receptor Tyrosine-based Inhibitory Motifs" (consensus ITIM: [I/V/L]-x-pY-x-x-[I/V/L]; death receptor variant: A-x-pY-x-x-L). This binding activates Shp-1 phosphatase by displacing an inhibitory interaction between the N-SH2 and PTP domains. By dephosphorylating the downstream elements of various signal transduction pathways, Shp-1 acts as a negative regulator. Shp-1 is highly expressed in hematopoietic cells and to a lesser extent in epithelial cells. It plays a significant role in various stages of hematopoietic development and in limiting neutrophil activation and inflammatory tissue damage through down-regulation of anti-apoptotic, cytokine-derived signals. In most lymphomas and leukemias Shp-1 expression is decreased or absent and Shp-1 may play a role in the pathogenesis of these and other cancers.

Citations: 2

[View Online »](#)

## Ordering Information

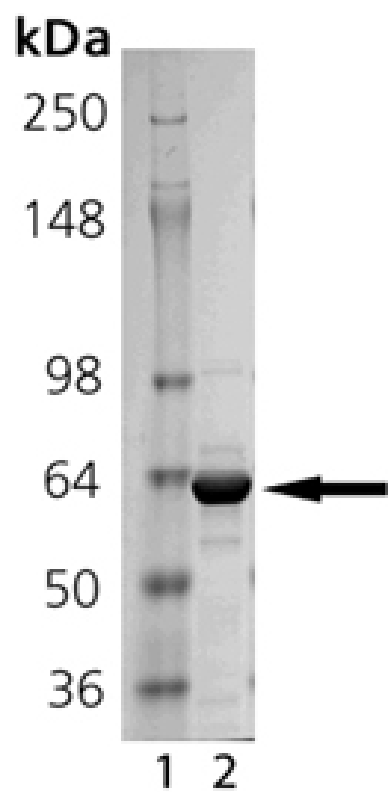
[Order Online »](#)

BML-SE334-0020

20µg

## Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW Marker; Lane 2: 2.0 $\mu$ g of purified Shp-1 (human) (recombinant).

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Alternative Name**      Src homology-2 containing protein tyrosine phosphatase 1

**Formulation**      Liquid. In 50mM TRIS/HCl, pH 8.0, containing 150mM NaCl, 5mM DTT, 0.03% Brij 35, 0.1mM EDTA and 10% glycerol.

**MW**      67.5 kDa

**Purity**      ≥90% (SDS-PAGE)

**Purity Detail**      Purified by multi-step chromatography.

**Source**      Produced in *E. coli*. Full length human Shp-1 (aa 1-595).

**Specific Activity**      ≥2000 pmol/min/μg assayed by p-nitrophenylphosphate (pNPP, 50 mM) hydrolysis at pH 7.0, 30°C

**UniProt ID**      P29350



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)