

# Shp-2 (human), (recombinant)

Shp-2 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. In most cases, full activation of the Erk/MAP kinase pathway by receptor tyrosine kinase signaling requires Shp-2, which is thought to act somewhere upstream of Ras.

## Ordering Information

[Order Online »](#)

<b>BML-SE273-0020</b>	20µg
-----------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	After initial defrosting, aliquot product into individual tubes and snap freeze, for example in liquid nitrogen or dry ice/ethanol.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Src homology-2 containing phosphatase 2
<b>Formulation</b>	Liquid. In 10mM MES, pH 5.7, containing 100mM NaCl, 5mM DTT and 33% glycerol.
<b>MW</b>	68 kDa
<b>Purity</b>	≥90% (SDS-PAGE)
<b>Purity Detail</b>	Purified by multi-step chromatography.
<b>Source</b>	Produced in <i>E. coli</i> . full length human Shp-2 (aa 1-593).
<b>Specific Activity</b>	≥100pmol/min/μg assayed by p-nitrophenylphosphate (pNPP, 50mM) hydrolysis at pH 7.0, 30°C.
<b>UniProt ID</b>	Q06124



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)