

# SHPS-1/SIRP-1 $\alpha$ polyclonal antibody

SHPS-1 is a member of the gene family called the signal regulatory proteins of which there are at least fifteen members. SHPS-1 is a substrate of many activated tyrosine kinases such as Insulin receptor and EGFR, amongst others. SHPS-1 has regulatory effects on cellular responses induced by serum, growth factors, insulin, oncogenes, growth hormones and cell adhesion and plays a general role in different physiological and pathological processes.

This antibody is covered by our [Worry-Free Guarantee](#).

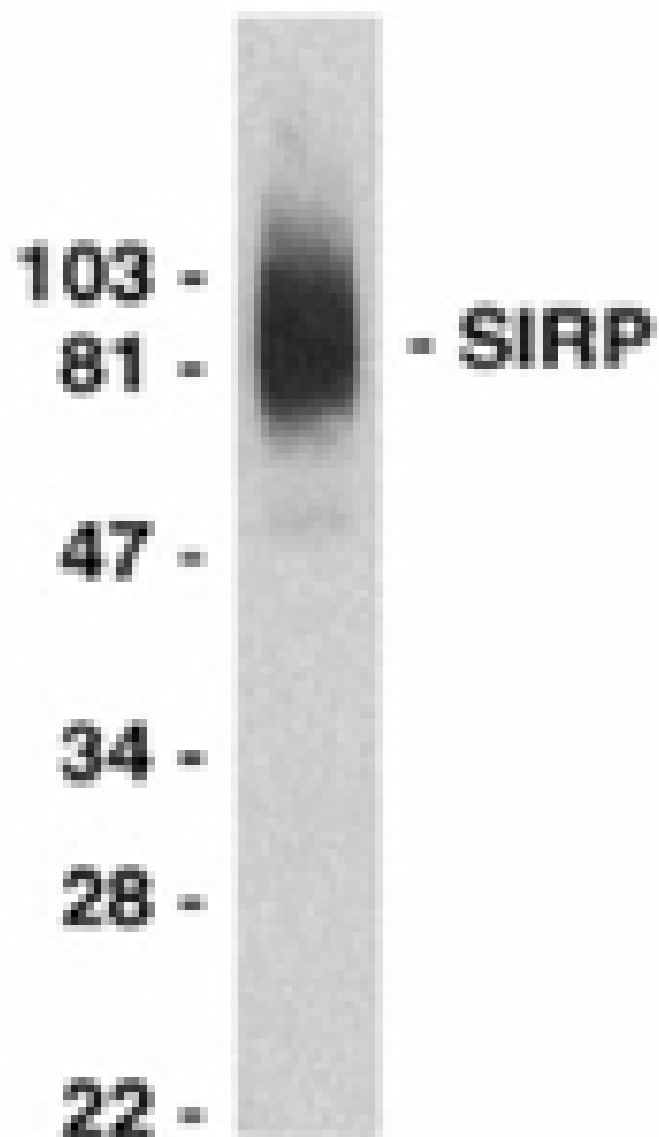
## Ordering Information

[Order Online »](#)

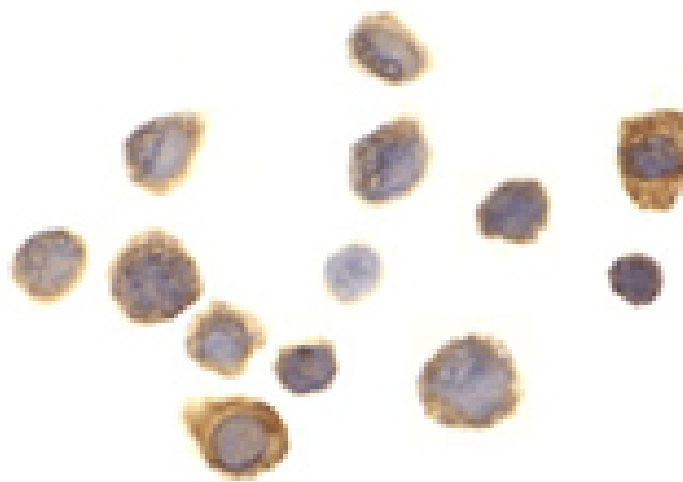
ADI-KAP-TK140-E	100 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of SIRP alpha in THP-1 whole cell lysate with SIRP alpha antibody at 1:2000 dilution.



Immunocytochemistry analysis of SIRP alpha in THP-1 cells with SIRP alpha antibody at 1 $\mu$ g/ml.

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application ICC, WB

Application Notes Detects a band of ~80-95kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

GenBank ID D86043

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human SHPS-1/SIRP-1 $\alpha$ .

Purity Detail Affinity chromatography purified via peptide column.

**Recommendation**  
**Dilutions/Conditions** Immunocytochemistry (1 $\mu$ g/ml)Western Blot (0.5-1.0 $\mu$ g/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID P78324

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



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)