SHPS-1/SIRP-1 α 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

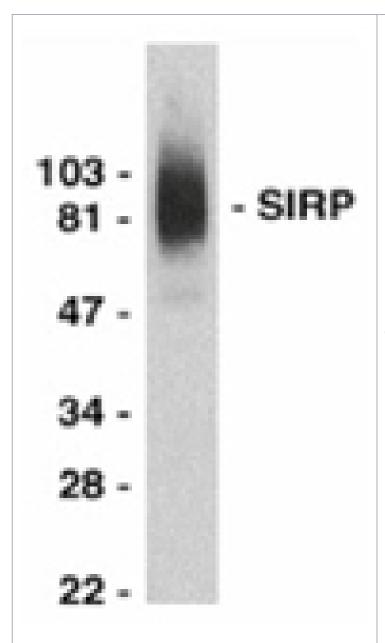
Order Online »

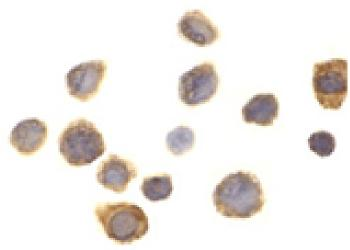
ADI-KAP-TK140-E

100µg

Manuals, SDS & CofA

View Online »





Immunocytochemistry analysis of SIRP alpha in THP-1 cells with SIRP alpha antibody at 1µg/ml.

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

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, WB

Application NotesDetects 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-1α.

Purity Detail Affinity chromatography purified via peptide column.

Recommendation Dilutions/Conditions Immunocytochemistry (1µg/ml)Western Blot (0.5-1.0µg/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

•

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com