

SHISA5 Polyclonal Antibody

(Catalog # A60898)

Background

Can induce apoptosis in a caspase-dependent manner and plays a role in p53/TP53-dependent apoptosis.

Description

SHISA5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q8N114

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein shisa-5 protein (56-137AA)

Storage

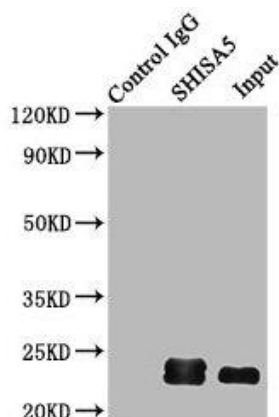
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

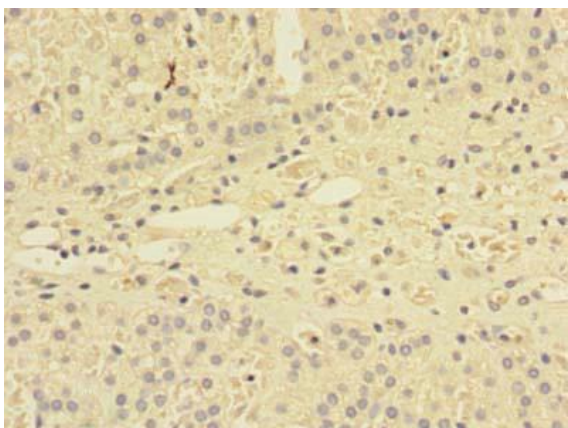
Protein shisa-5, SHISA5, SCOTIN, PSEC0133, Putative NF-kappa-B-activating protein 120, Scotin

Application

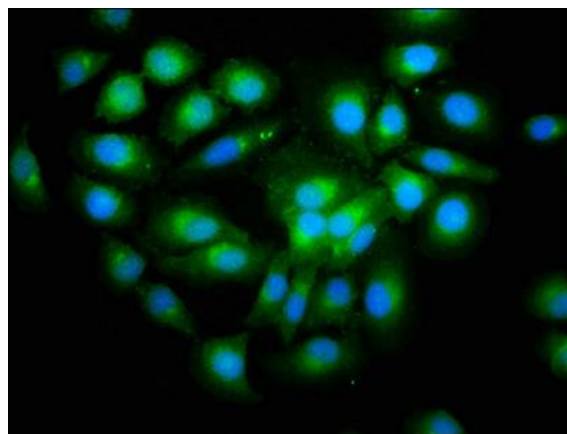
ELISA, IHC, IF, IP; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:500, IP:1:200-1:2000



Immunoprecipitating SHISA5 in HepG2 whole cell lysate
Lane 1: Rabbit control IgG (1ug) instead of SHISA5 Ab in HepG2 whole cell lysate.
For western blotting, an HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000)
Lane 2: SHISA5 Ab (6ug) + HepG2 whole cell lysate (500ug)
Lane 3: HepG2 whole cell lysate (10ug)



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using SHISA5 Antibody at dilution of 1:100



Immunofluorescence staining of A549 cells with SHISA5 Antibody at 1:200, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).