

SIPA1L2 Polyclonal Antibody

(Catalog #A60914)

Description

SIPA1L2 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

Q9P2F8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Signal-induced proliferation-associated 1-like protein 2 protein (522-778AA)

Storage

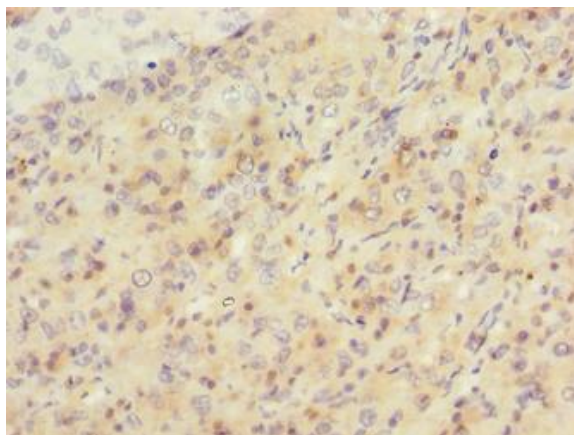
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

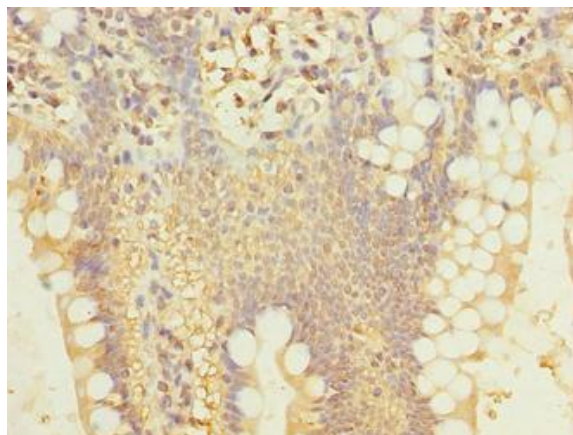
Signal-induced proliferation-associated 1-like protein 2, SIPA1-like protein 2, SIPA1L2, KIAA1389

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human liver cancer using SIPA1L2 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human small intestine tissue using SIPA1L2 Polyclonal Antibody at dilution of 1:100