
FARP1 Polyclonal Antibody

(Catalog # A68226)

Background

Functions as guanine nucleotide exchange factor for RAC1. May play a role in semaphorin signaling. Plays a role in the assembly and disassembly of dendritic filopodia, the formation of dendritic spines, regulation of dendrite length and ultimately the formation of synapses.

Description

FARP1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9Y4F1

Purification

Protein G purified

Immunogen

Recombinant Human FERM, ARHGEF and pleckstrin domain-containing protein 1 protein (1-300AA)

Storage

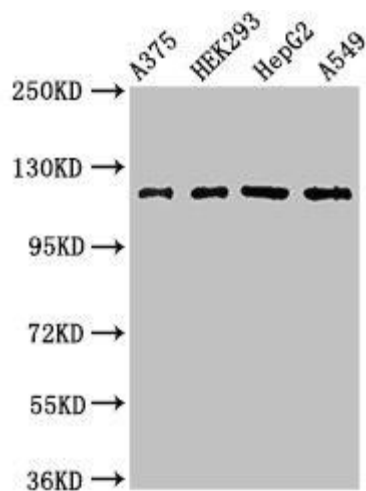
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

FERM, RhoGEF and pleckstrin domain-containing protein 1, Chondrocyte-derived ezrin-like protein, Pleckstrin homology domain-containing family C member 2, PH domain-containing family C member 2, FARP1, CDEP, PLEKHC2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: A375 whole cell lysate, HEK293 whole cell lysate, HepG2 whole cell lysate, A549 whole cell lysate

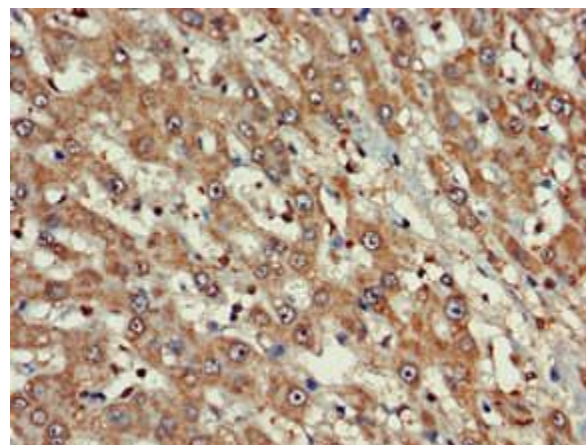
All lanes: FARP1 Antibody at 3µg/ml

Secondary

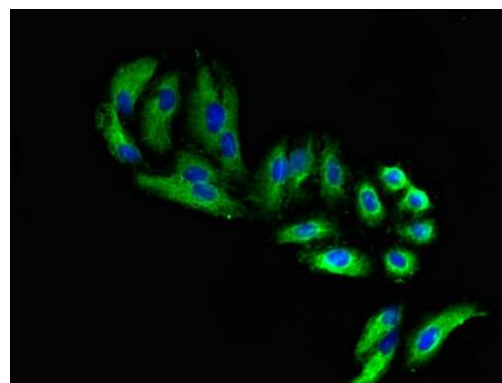
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 119, 123, 15 kDa

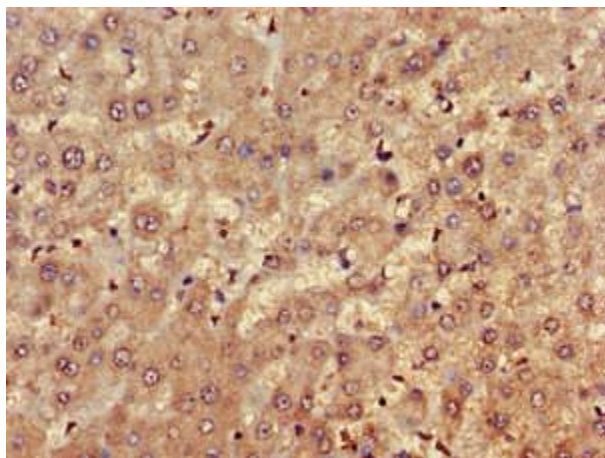
Observed band size: 119 kDa



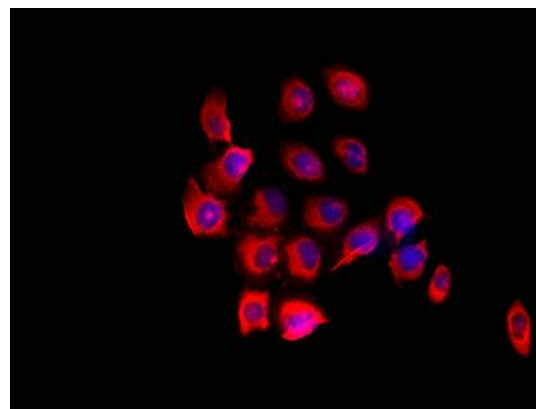
Immunohistochemistry of paraffin-embedded human liver cancer using FARP1 Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using FARP1 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human liver tissue using FARP1 Antibody at dilution of 1:100



Immunofluorescent analysis of A549 cells using FARP1 Antibody at dilution of 1:100 and Cy3-conjugated Goat Anti-Rabbit IgG