
RAP1A Polyclonal Antibody

(Catalog # A60518)

Background

Induces morphological reversion of a cell line transformed by a Ras oncogene. Counteracts the mitogenic function of Ras, at least partly because it can interact with Ras GAPs and RAF in a competitive manner. Together with ITGB1BP1, regulates KRIT1 localization to microtubules and membranes. Plays a role in nerve growth factor (NGF)-induced neurite outgrowth. Plays a role in the regulation of embryonic blood vessel formation. Involved in the establishment of basal endothelial barrier function. May be involved in the regulation of the vascular endothelial growth factor receptor KDR expression at endothelial cell-cell junctions.

Description

RAP1A 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

P62834

Purification

Protein G purified

Immunogen

Recombinant Human Ras-related protein Rap-1A protein (1-181AA)

Storage

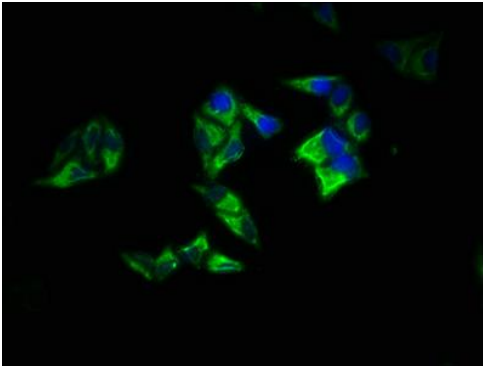
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

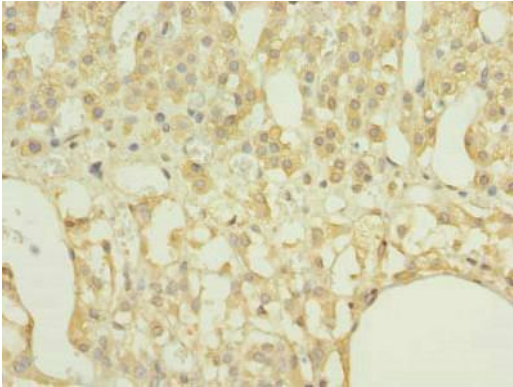
Ras-related protein Rap-1A, RAP1A, KREV1, C21KG, G-22K, GTP-binding protein smg p21A, Ras-related protein Krev-1

Application

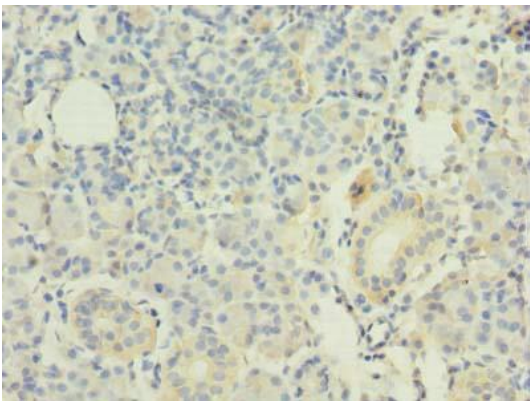
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:500



Immunofluorescence staining of HeLa cells with RAP1A 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).



Immunohistochemistry of paraffin-embedded human pancreatic tissue using RAP1A Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human pancreatic tissue using RAP1A Antibody at dilution of 1:100