
PIK3AP1 Polyclonal Antibody

(Catalog #A68967)

Background

Signaling adapter that contributes to B-cell development by linking B-cell receptor (BCR) signaling to the phosphoinositide 3-kinase (PI3K)-Akt signaling pathway. Has a complementary role to the BCR coreceptor CD19, coupling BCR and PI3K activation by providing a docking site for the PI3K subunit PIK3R1. Alternatively, links Toll-like receptor (TLR) signaling to PI3K activation, a process preventing excessive inflammatory cytokine production. Also involved in the activation of PI3K in natural killer cells. May be involved in the survival of mature B-cells via activation of REL.

Description

PIK3AP1 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

Q6ZUJ8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Phosphoinositide 3-kinase adapter protein 1 protein (293-404AA)

Storage

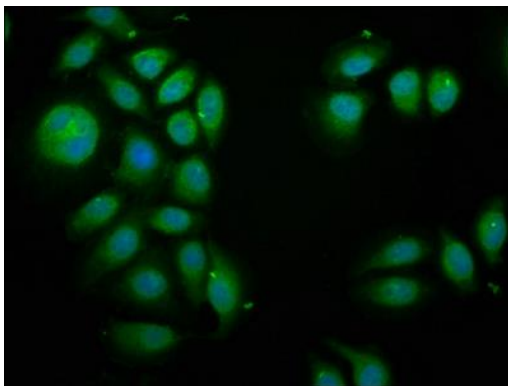
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Phosphoinositide 3-kinase adapter protein 1, B-cell adapter for phosphoinositide 3-kinase, B-cell phosphoinositide 3-kinase adapter protein 1, PIK3AP1, BCAP

Application

ELISA, IF; Recommended dilution: IF 1:50 - 1:200



Immunofluorescence staining of A549 cells with PIK3AP1 Polyclonal Antibody at 1:100, 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).