
PLPP3 Polyclonal Antibody

(Catalog # A63190)

Background

Catalyzes the conversion of phosphatidic acid (PA) to diacylglycerol (DG). In addition, it hydrolyzes lysophosphatidic acid (LPA), ceramide-1-phosphate (C-1-P) and sphingosine-1-phosphate (S-1-P). The relative catalytic efficiency is LPA = PA > C-1-P > S-1-P. May be involved in cell adhesion and in cell-cell interactions.

Description

PLPP3 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

O14495

Purification

Protein G purified

Immunogen

Recombinant Human Phospholipid phosphatase 3 protein (144-193AA)

Storage

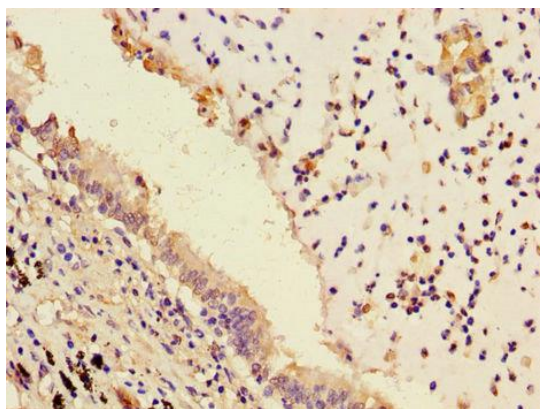
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

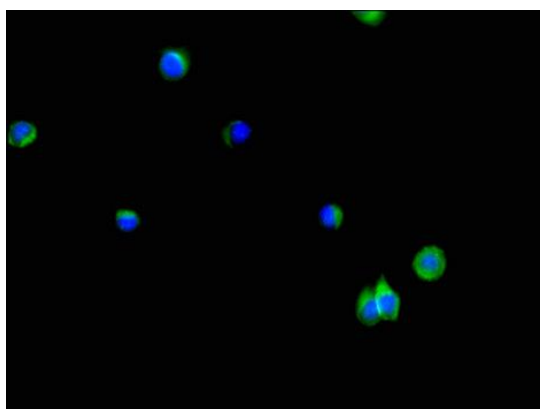
Lipid phosphate phosphohydrolase 3, PAP2-beta, Phosphatidate phosphohydrolase type 2b, Phosphatidic acid phosphatase 2b, PAP-2b, PAP2b, Vascular endothelial growth factor and type I collagen-inducible protein, VCIP, PPAP2B, LPP3

Application

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



Immunohistochemistry of paraffin-embedded human lung cancer tissue at dilution of 1:100



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