
PPAP2B Polyclonal Antibody

(Catalog #A72517)

Background

The protein encoded by this gene is a member of the phosphatidic acid phosphatase (PAP) family. PAPs convert phosphatidic acid to diacylglycerol, and function in de novo synthesis of glycerolipids as well as in receptor-activated signal transduction mediated by phospholipase D. This protein is a membrane glycoprotein localized at the cell plasma membrane. It has been shown to actively hydrolyze extracellular lysophosphatidic acid and short-chain phosphatidic acid. The expression of this gene is found to be enhanced by epidermal growth factor in Hela cells.

Description

PPAP2B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

O14495

Purification

Affinity Purification

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 1-100 of human PPAP2B (NP_003704.3).

Storage

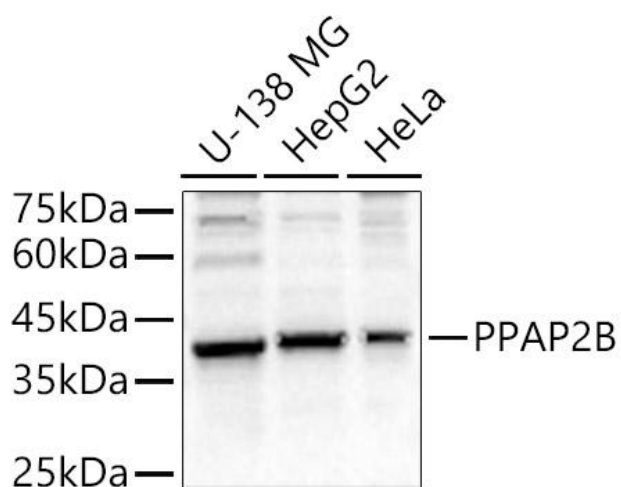
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

PLPP3; Dri42; LPP3; PAP2B; PPAP2B; VCIP; phospholipid phosphatase 3

Application

WB, IHC; Recommended dilution: WB 1:500 - 1:1000, IHC 1:100 - 1:200



Western blot analysis of various lysates, using PPAP2B Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25µg per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 10s.