

PLA2G7 Polyclonal Antibody

(Catalog # A55383)

Background

Modulates the action of platelet-activating factor (PAF) by hydrolyzing the sn-2 ester bond to yield the biologically inactive lyso-PAF. Has a specificity for substrates with a short residue at the sn-2 position. It is inactive against long-chain phospholipids.

Description

PLA2G7 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q13093

Purification

Protein G purified

Immunogen

Recombinant Human Platelet-activating factor acetylhydrolase protein (49-441AA)

Storage

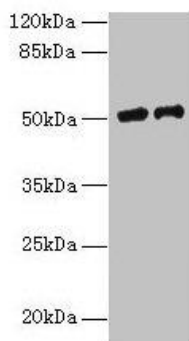
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

PAF acetylhydrolase, 1-alkyl-2-acetyl-glycerophosphocholine esterase, 2-acetyl-1-alkyl-glycerophosphocholine esterase, Group-VIIA phospholipase A2, gVIIA-PLA2, LDL-associated phospholipase A2, LDL-PLA(2), PAF 2-acylhydrolase, PLA2G7, PAFAH

Application

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



Western blot

All lanes: PLA2G7 antibody at 9ug/ml

Lane 1: RAW264.7 whole cell lysate

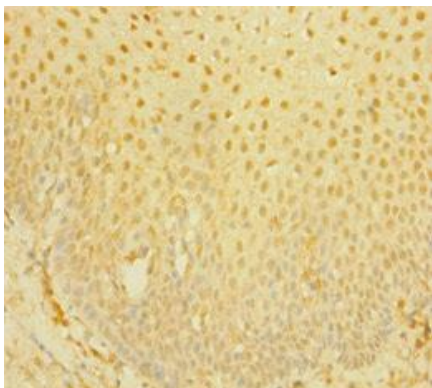
Lane 2: Mouse liver tissue

Secondary

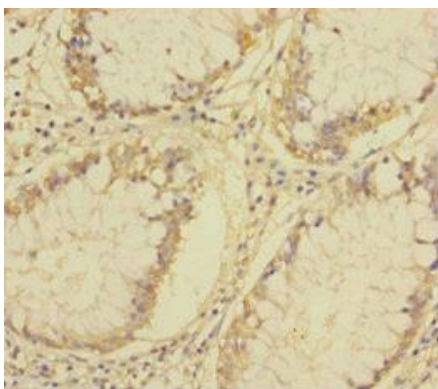
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 51 kDa

Observed band size: 51 kDa



Immunohistochemistry of paraffin-embedded human tonsil tissue using PLA2G7 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human colon cancer using PLA2G7 Antibody at dilution of 1:100