
PLA1A Polyclonal Antibody

(Catalog # A69655)

Background

Hydrolyzes the ester bond at the sn-1 position of glycerophospholipids and produces 2-acyl lysophospholipids. Hydrolyzes phosphatidylserine (PS) in the form of liposomes and 1-acyl-2 lysophosphatidylserine (lyso-PS), but not triolein, phosphatidylcholine (PC), phosphatidylethanolamine (PE), phosphatidic acid (PA) or phosphatidylinositol (PI). Isoform 2 hydrolyzes lyso-PS but not PS. Hydrolysis of lyso-PS in peritoneal mast cells activated by receptors for IgE leads to stimulate histamine production.

Description

PLA1A 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

Q53H76

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Phospholipase A1 member A protein (293-398AA)

Storage

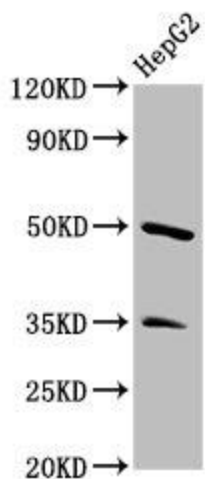
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Phospholipase A1 member A, Phosphatidylserine-specific phospholipase A1, PS-PLA1, PLA1A, NMD, PSPLA1

Application

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



Western Blot

Positive WB detected in: HepG2 whole cell lysate

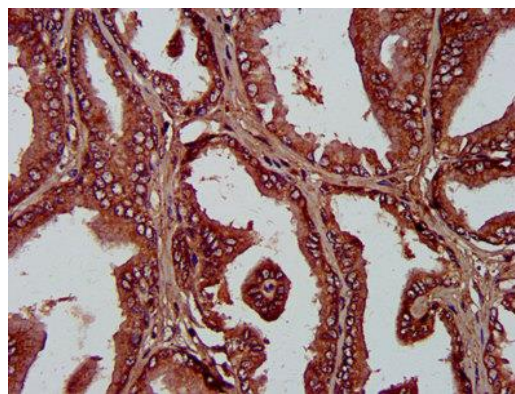
All lanes: PLA1A Polyclonal Antibody at 3.4ug/ml

Secondary

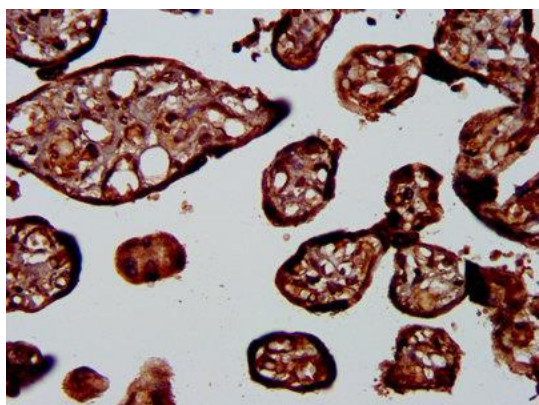
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 50, 41, 48, 32 KDa

Observed band size: 50 KDa



IHC image of PLA1A Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human prostate tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of PLA1A Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human placenta tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.\