

PNPLA1 Polyclonal Antibody

(Catalog # A66754)

Background

Lipid hydrolase. Important in the formation of the epidermal lipid barrier. Plays a role in glycerophospholipid metabolism.

Description

PNPLA1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q8N8W4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Patatin-like phospholipase domain-containing protein 1 protein (101-300AA)

Storage

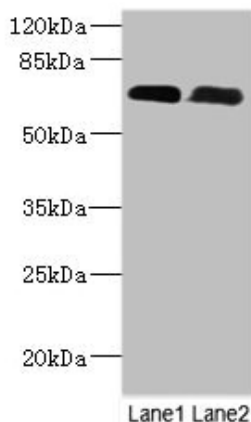
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Patatin-like phospholipase domain-containing protein 1, PNPLA1

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

All lanes: PNPLA1 Polyclonal Antibody at 12ug/ml

Lane 1: Mouse spleen tissue

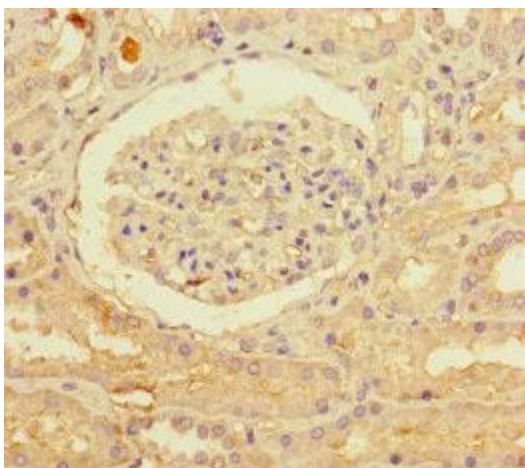
Lane 2: Mouse skeletal muscle tissue

Secondary

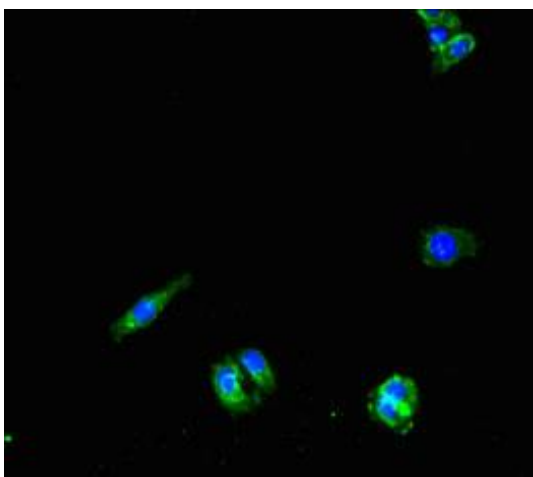
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 58, 48, 49 kDa

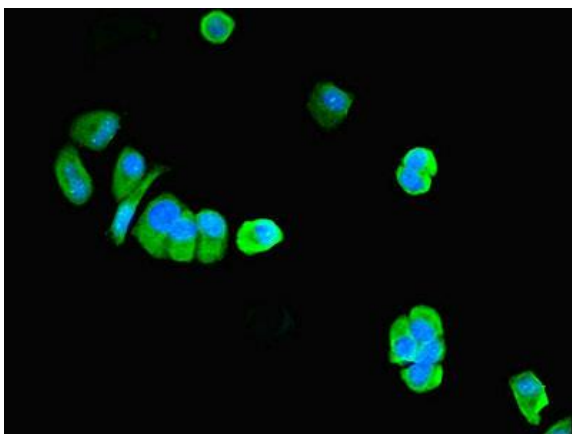
Observed band size: 58 kDa



Immunohistochemistry of paraffin-embedded human kidney tissue using PNPLA1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG-2 cells using PNPLA1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunofluorescent analysis of MCF-7 cells using PNPLA1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)