

PLA2G2A Polyclonal Antibody

(Catalog # A52245)

Background

Thought to participate in the regulation of the phospholipid metabolism in biomembranes including eicosanoid biosynthesis. Catalyzes the calcium-dependent hydrolysis of the 2-acyl groups in 3-sn-phosphoglycerides.

Description

PLA2G2A 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

P14555

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Phospholipase A2, membrane associated protein (21-144AA)

Storage

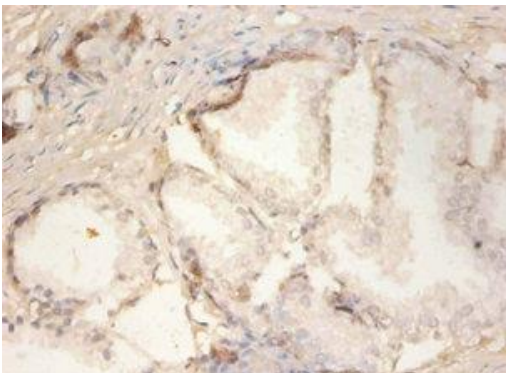
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

GIIC sPLA2, Group IIA phospholipase A2, Non-pancreatic secretory phospholipase A2, Phosphatidylcholine 2-acylhydrolase 2A, PLA2G2A, PLA2B, PLA2L, RASF-A

Application

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



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