
SMPD3 Polyclonal Antibody

(Catalog #A67638)

Background

Catalyzes the hydrolysis of sphingomyelin to form ceramide and phosphocholine. Ceramide mediates numerous cellular functions, such as apoptosis and growth arrest, and is capable of regulating these 2 cellular events independently. Also hydrolyzes sphingosylphosphocholine. Regulates the cell cycle by acting as a growth suppressor in confluent cells. Probably acts as a regulator of postnatal development and participates in bone and dentin mineralization.

Description

SMPD3 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

Q9NY59

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Sphingomyelin phosphodiesterase 3 protein (401-655AA)

Storage

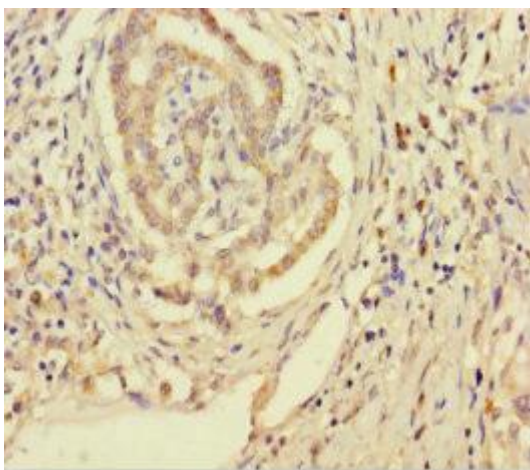
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

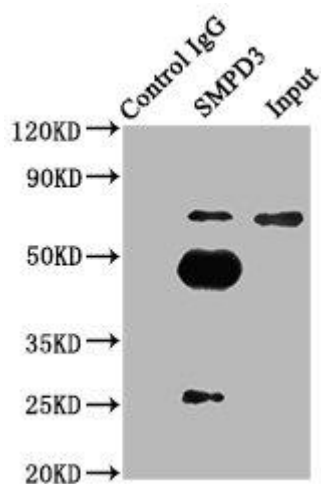
Sphingomyelin phosphodiesterase 3, Neutral sphingomyelinase 2, nSMase-2, nSMase2, Neutral sphingomyelinase II, SMPD3

Application

ELISA, WB, IHC, IP; Recommended dilution: WB: 1:500-1:2000, IHC:1:50-1:200, IP:1:200-1:2000



Immunohistochemistry of paraffin-embedded human pancreatic cancer using SMPD3 Antibody at dilution of 1:100



Immunoprecipitating SMPD3 in mouse brain whole cell lysate
 Lane 1: Rabbit control IgG (1 μ g) instead of SMPD3 Antibody in mouse brain whole cell lysate. For western blotting, a HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000)
 Lane 2: SMPD3 Antibody (6 μ g) + Mouse brain whole cell lysate (500 μ g)
 Lane 3: Mouse brain whole cell lysate (10 μ g)