

ANO4 Polyclonal Antibody

(Catalog # A64886)

Background

Has calcium-dependent phospholipid scramblase activity; scrambles phosphatidylserine, phosphatidylcholine and galactosylceramide. Does not exhibit calcium-activated chloride channel (CaCC) activity.

Description

ANO4 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

Q32M45

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Anoctamin-4 protein (1-200AA)

Storage

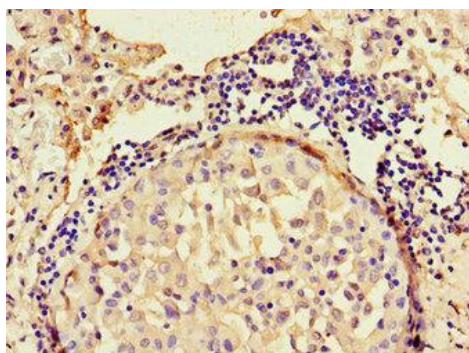
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

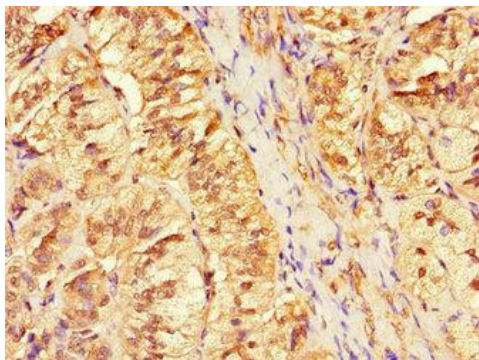
Anoctamin-4, Transmembrane protein 16D, ANO4, TMEM16D

Application

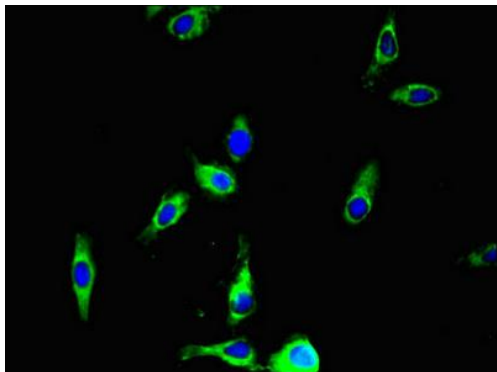
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human lung cancer using ANO4 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using ANO4 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using ANO4 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)