

PKD1L2 Polyclonal Antibody

(Catalog # A60286)

Background

May function as an ion-channel regulator. May function as a G-protein-coupled receptor.

Description

PKD1L2 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

Q7Z442

Purification

Affinity purified

Immunogen

Recombinant Human Polycystic kidney disease protein 1-like 2 protein (1-306AA)

Storage

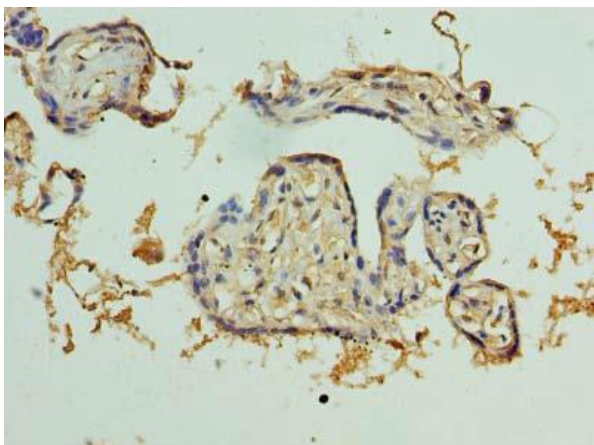
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

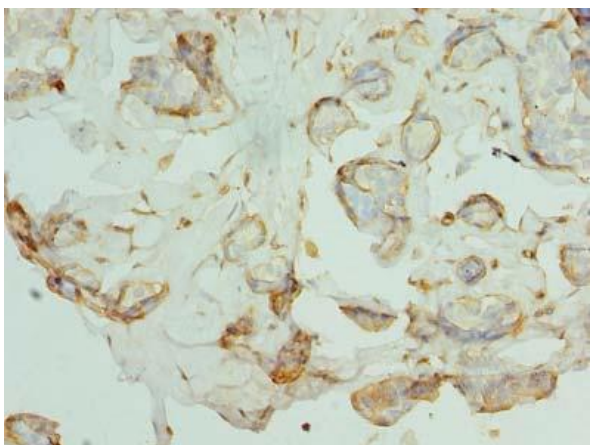
Polycystic kidney disease protein 1-like 2, PKD1L2, KIAA1879, PC1L2, PC1-like 2 protein, Polycystin-1L2

Application

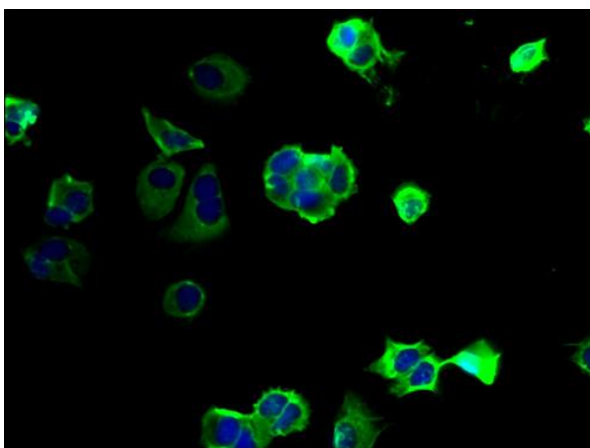
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human placenta tissue using PKD1L2 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human breast cancer using PKD1L2 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells using PKD1L2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)