

TMEM95 Polyclonal Antibody

(Catalog # A63676)

Description

TMEM95 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

Q3KNT9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transmembrane protein 95 protein (17-145AA)

Storage

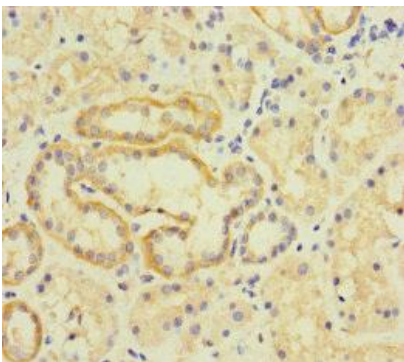
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

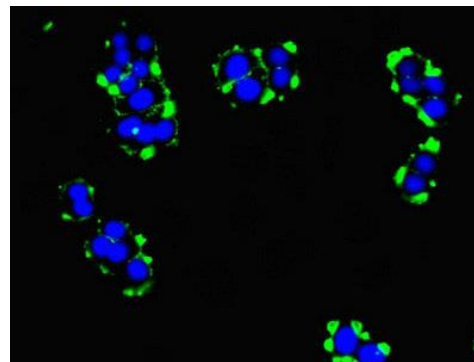
Transmembrane protein 95, TMEM95, UNQ9390/PRO34281

Application

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



Immunohistochemistry of paraffin-embedded human kidney tissue using TMEM95 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of PC3 cells using TMEM95 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)