

TMEM63A Polyclonal Antibody

(Catalog # A61442)

Background

Acts as an osmosensitive calcium-permeable cation channel.

Description

TMEM63A 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

O94886

Purification

>95%, Protein G purified

Immunogen

Recombinant Human CSC1-like protein 1 protein (212-420AA)

Storage

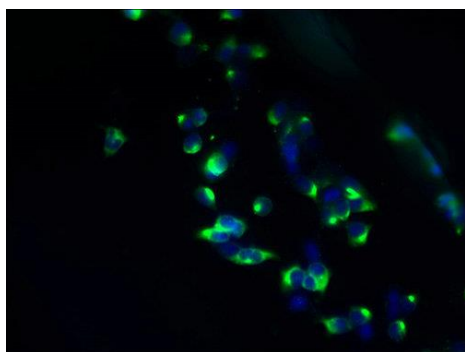
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

CSC1-like protein 1, TMEM63A, KIAA0489, KIAA0792, Transmembrane protein 63A

Application

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



Immunofluorescent analysis of 293T cells at a dilution of 1:100
and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit
IgG(H+L)