

TMEM47 Polyclonal Antibody

(Catalog # A63644)

Background

cell junction, plasma membrane, cell-cell adhesion

Description

TMEM47 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

Q9BQJ4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transmembrane protein 47 protein (42-82AA)

Storage

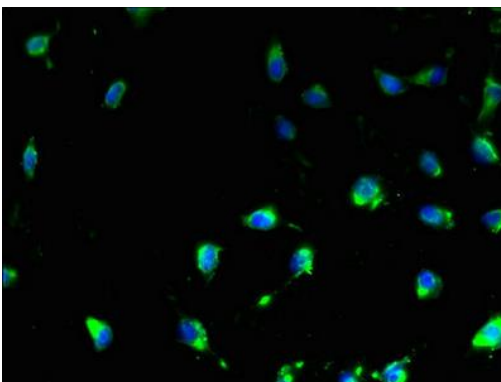
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

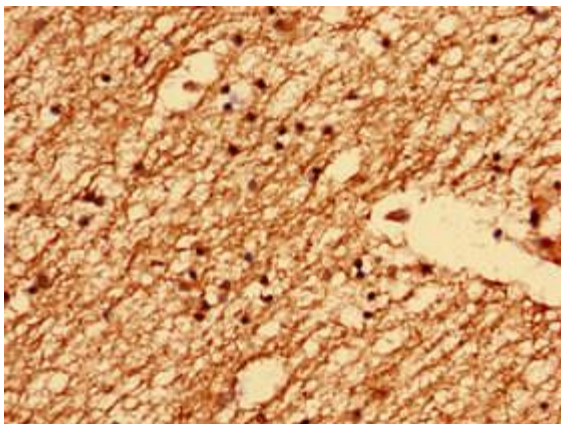
Transmembrane protein 47, Brain cell membrane protein 1, Transmembrane 4 superfamily member 10, TMEM47, BCMP1, TM4SF10

Application

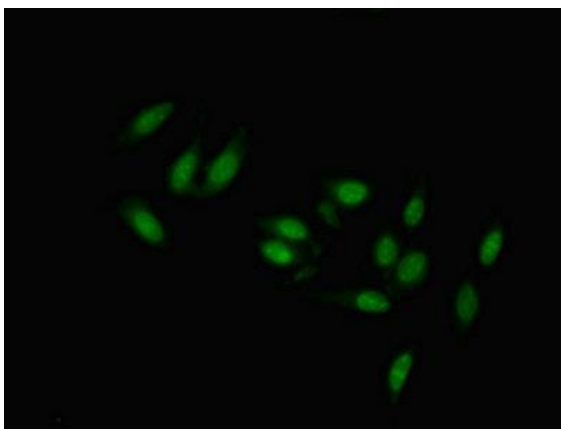
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunofluorescent analysis of U251 cells using TMEM47 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunohistochemistry analysis of human brain tissue using TMEM47 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using TMEM47 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)