

TSPAN14 Polyclonal Antibody

(Catalog # A61554)

Background

integral component of plasma membrane, plasma membrane, enzyme binding, establishment of protein localization to plasma membrane, positive regulation of Notch signaling pathway, protein maturation

Description

TSPAN14 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

Q8NG11

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tetraspanin-14 protein (114-215AA)

Storage

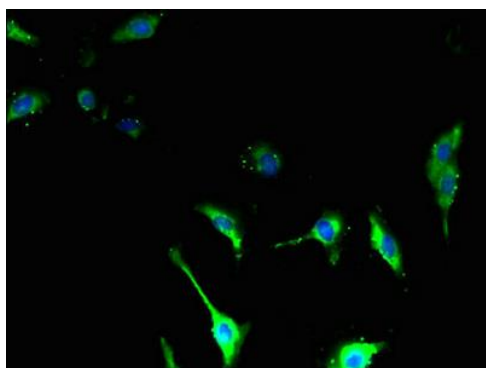
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

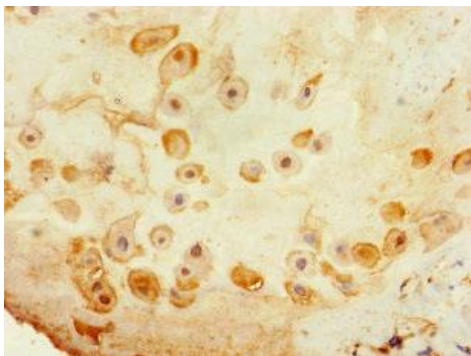
Tetraspanin-14, Tspan-14, TSPAN14, TM4SF14, DC-TM4F2, Transmembrane 4 superfamily member 14

Application

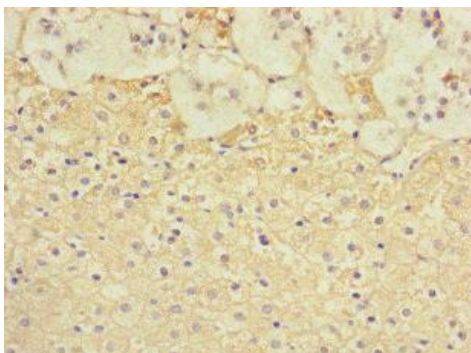
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunofluorescent analysis of U251 cells using TSPAN14 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human placenta tissue using TSPAN14 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using TSPAN14 Antibody at dilution of 1:100