

TSPAN16 Polyclonal Antibody

(Catalog # A61562)

Background

Integral component of membrane, integral component of plasma membrane, cell surface receptor signaling pathway

Description

TSPAN16 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

Q9UKR8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tetraspanin-16 protein (116-160AA)

Storage

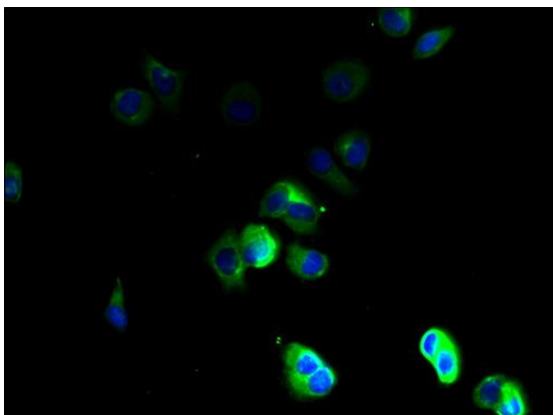
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Tetraspanin-16, Tspan-16, TSPAN16, TM4SF16, Tetraspanin TM4-B, Transmembrane 4 superfamily member 16

Application

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



Immunofluorescent analysis of MCF-7 cells using TSPAN16 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)