

TSPAN5 Polyclonal Antibody

(Catalog # A61566)

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

TSPAN5 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

P62079

Purification

Affinity purified

Immunogen

Recombinant Human Tetraspanin-5 protein (114-232AA)

Storage

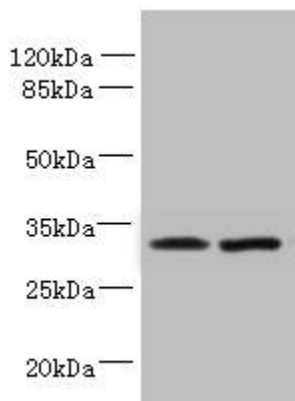
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Tetraspanin-5, Tspan-5, TSPAN5, TM4SF9, Tetraspan NET-4, Transmembrane 4 superfamily member 9

Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200



Western blot

All lanes: TSPAN5 Polyclonal Antibody at 1.2ug/ml

Lane 1: Jurkat whole cell lysate

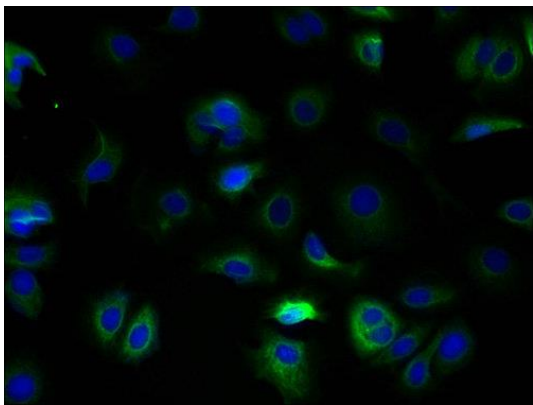
Lane 2: A549 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 30 kDa

Observed band size: 30 kDa



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