

TSPAN31 Polyclonal Antibody

(Catalog # A50933)

Background

TSPAN31 (Tetraspanin 31) is a protein-coding gene that belongs to the transmembrane 4 superfamily, or tetraspanin family of proteins. Most of these members are cell-surface proteins and are characterized by the presence of four hydrophobic domains. They are known to mediate signal transduction events that play a role in the regulation of cell development, activation, growth, and motility. This encoded protein is thought to be involved in growth-related cellular processes. TSPAN31 is associated with tumorigenesis and osteosarcoma. Diseases associated with this gene include Well-Differentiated Liposarcoma and Familial Retinoblastoma.

Description

TSPAN31 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

Q12999

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tetraspanin-31 protein (94-173AA)

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

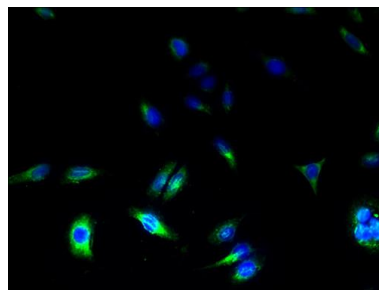
Tetraspanin-31, Tspan-31, Sarcoma-amplified sequence, TSPAN31, SAS

Application

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



Immunohistochemistry of paraffin-embedded human placenta tissue using TSPAN31 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using TSPAN31 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)