

TMEM53 Polyclonal Antibody

(Catalog # A63648)

Background

Located in the nucleus. Ensures normal bone formation through the negative regulation of bone morphogenetic protein (BMP) signaling in osteoblast lineage cells by blocking cytoplasm-nucleus translocation of phosphorylated SMAD1/5/9 proteins.

Description

TMEM53 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q6P2H8

Purification

Protein G purified

Immunogen

Recombinant Human Transmembrane protein 53 protein (1-170AA)

Storage

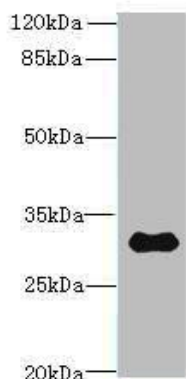
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transmembrane protein 53, TMEM53

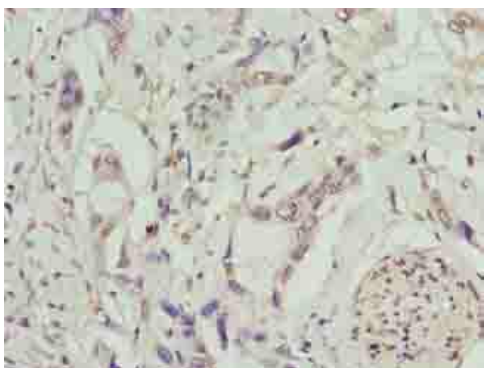
Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200

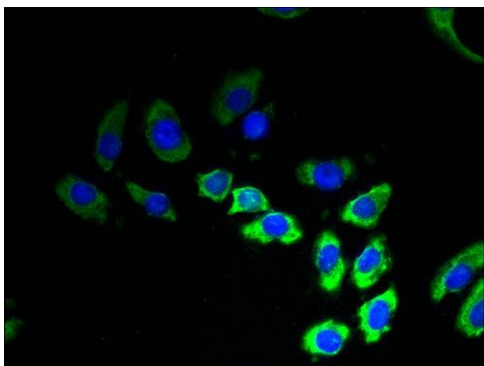


Western blot

All lanes: TMEM53 Polyclonal Antibody at
5ug/ml+Mouse brain tissue
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 32, 24 kDa
Observed band size: 32 kDa



Immunohistochemistry of paraffin-embedded human pancreatic cancer at dilution of 1:100



Immunofluorescent analysis of A549 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)