

TMBIM1 Polyclonal Antibody

(Catalog # A61338)

Background

Negatively regulates aortic matrix metalloproteinase-9 (MMP9) production and may play a protective role in vascular remodeling.

Description

TMBIM1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q969X1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein lifeguard 3 protein (1-109AA)

Storage

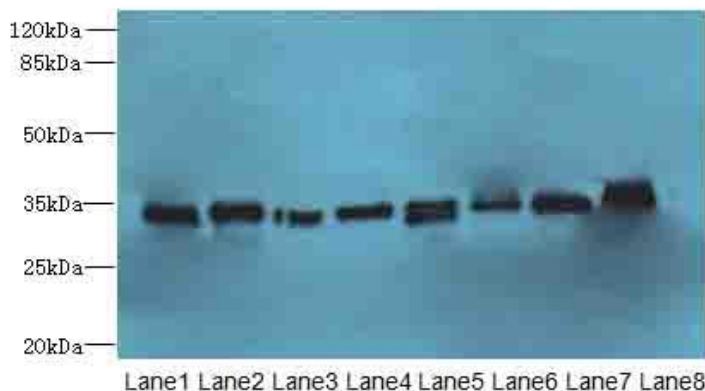
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein lifeguard 3, TMBIM1, LFG3, RECS1, PP1201, PSEC0158, Protein RECS1 homolog, Transmembrane BAX inhibitor motif-containing protein 1

Application

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



Western blot

All Lanes: TMBIM1 Polyclonal Antibody at 8ug/ml

Lane 1: U251 whole cell lysate

Lane 2: K562 whole cell lysate

Lane 3: A549 whole cell lysate

Lane 4: MCF7 whole cell lysate

Lane 5: Mouse lung tissue

Lane 6: Mouse liver tissue

Lane 7: Mouse kidney tissue

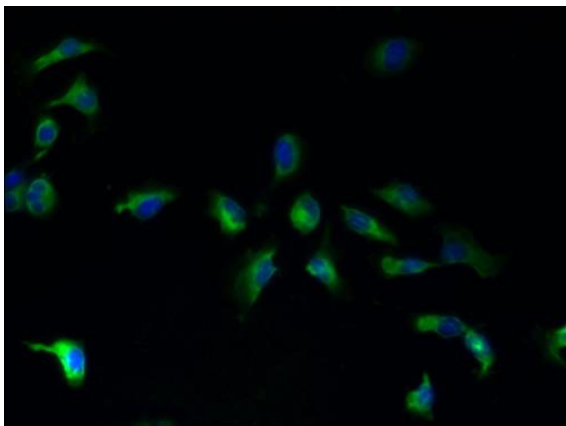
Lane 8: Human placenta tissue

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa



Immunofluorescent analysis of U251 cells using TMBIM1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)