

TMEM65 Polyclonal Antibody

(Catalog # A63656)

Description

TMEM65 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

Q6PI78

Purification

Affinity Purified

Immunogen

Recombinant Human Transmembrane protein 65 protein (63-240AA)

Storage

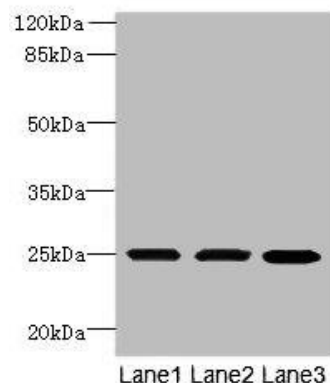
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

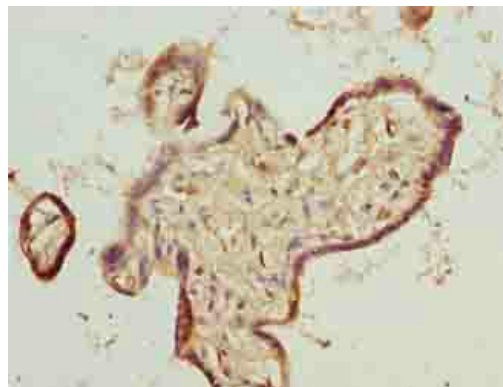
Transmembrane protein 65, TMEM65

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: TMEM65 Polyclonal Antibody at 6ug/ml
Lane 1: Mouse liver tissue, Lane 2: Mouse kidney tissue
Lane 3: Mouse gonadal tissue
Secondary: Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 25 kDa
Observed band size: 25 kDa



Immunohistochemistry of paraffin-embedded human placenta tissue at dilution of 1:100