

TMEM9 Polyclonal Antibody

(Catalog # A61446)

Background

May be involved in intracellular transport.

Description

TMEM9 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

Q9P0T7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transmembrane protein 9 protein (111-183AA)

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

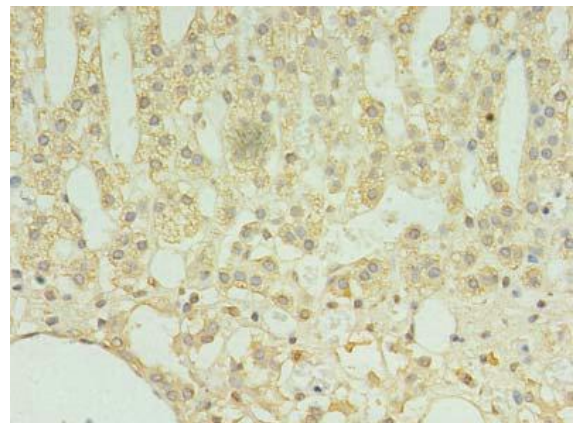
Transmembrane protein 9, TMEM9, DERP4, HSPC186, PSEC0012, UNQ631/PRO1248, Dermal papilla-derived protein 4

Application

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



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



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