

TMEM140 Polyclonal Antibody

(Catalog # A63584)

Description

TMEM140 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

Q9NV12

Purification

Affinity purified

Immunogen

Recombinant Human Transmembrane protein 140 protein (33-77AA)

Storage

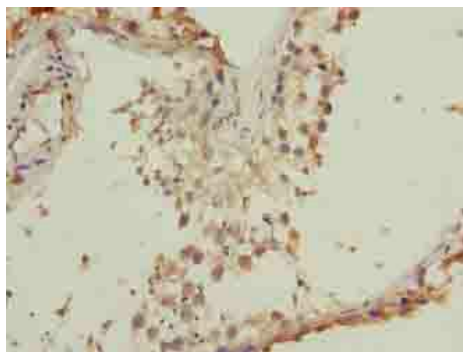
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

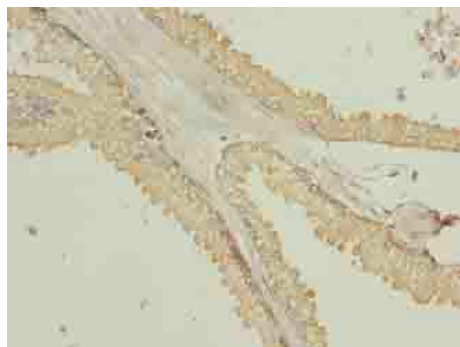
Transmembrane protein 140, TMEM140

Application

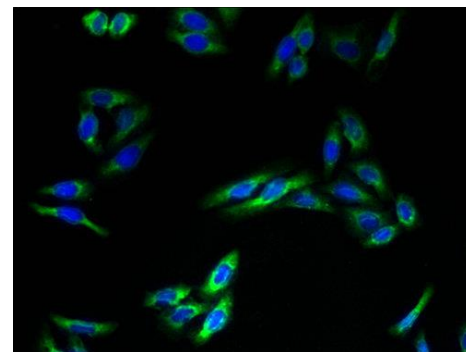
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



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



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



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