

## TM7SF3 Polyclonal Antibody

(Catalog # A61334)

### Background

extracellular exosome

### Description

TM7SF3 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

Q9NS93

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Transmembrane 7 superfamily member 3 protein (22-295AA)

### Storage

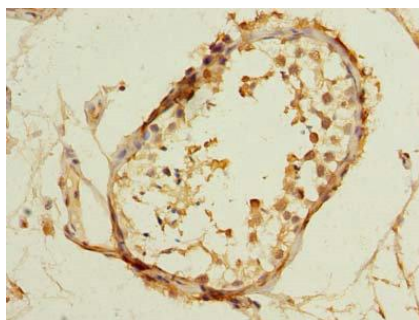
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

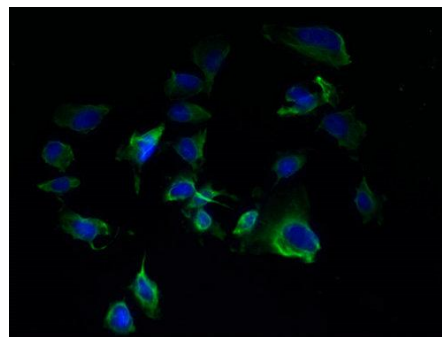
Transmembrane 7 superfamily member 3, TM7SF3, Seven span transmembrane protein

### Application

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



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



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