
TTC16 Polyclonal Antibody

(Catalog A61582)

Description

TTC16 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

Q8NEE8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tetratricopeptide repeat protein 16 protein (551-873AA)

Storage

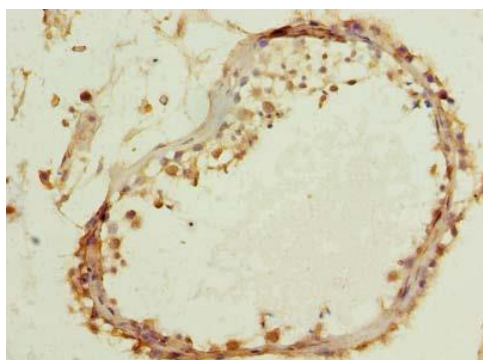
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Tetratricopeptide repeat protein 16, TPR repeat protein 16, TTC16

Application

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



Immunohistochemistry of paraffin-embedded human testis tissue using TTC16 Polyclonal Antibody at dilution of 1:100