

## TMTC4 Polyclonal Antibody

(Catalog # A61454)

### Background

Transfers mannosyl residues to the hydroxyl group of serine or threonine residues. The 4 members of the TMTC family are O-mannosyl-transferases dedicated primarily to the cadherin superfamily; each member seems to have a distinct role in decorating the cadherin domains with O-linked mannose glycans at specific regions. Also acts as O-mannosyl-transferase on other proteins such as PDIA3.

### Description

TMTC4 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

Q5T4D3

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Transmembrane and TPR repeat-containing protein 4 protein (245-376AA)

### Storage

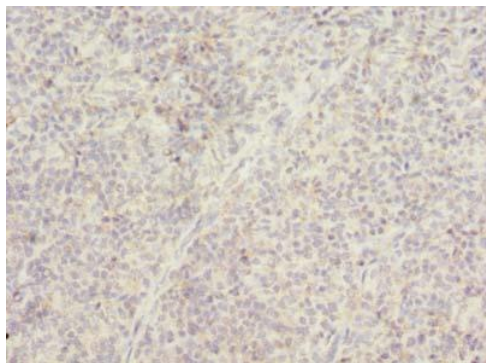
Shipped at 4°C. Upon delivery, aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Transmembrane and TPR repeat-containing protein 4, TMTC4

### Application

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



Immunohistochemistry of paraffin-embedded human tonsil using TMTC4 Polyclonal Antibody at dilution of 1:100 the input