

## TTC5 Polyclonal Antibody

(Catalog # A66330)

### Background

Adapter protein involved in p53/TP53 response that acts by regulating and mediating the assembly of multi-protein complexes. Required to facilitate the interaction between JMY and p300/EP300 and increase p53/TP53-dependent transcription and apoptosis. Prevents p53/TP53 degradation by MDM2.

### Description

TTC5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

Q8N0Z6

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Tetratricopeptide repeat protein 5 protein (201-440AA)

### Storage

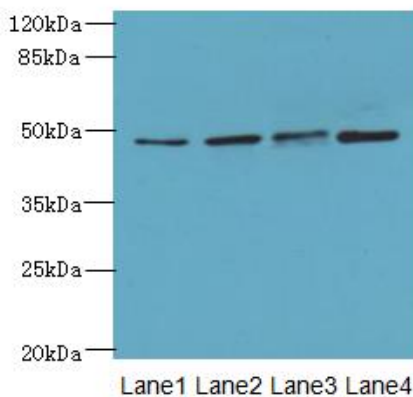
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Tetratricopeptide repeat protein 5, TPR repeat protein 5, Stress-responsive activator of p300, Strap, TTC5

### Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



### Western blot

All Lanes: TTC5 Polyclonal Antibody at 6ug/ml

Lane 1: HepG-2 whole cell lysate

Lane 2: A549 whole cell lysate

Lane 3: MCF7 whole cell lysate

Lane 4: Mouse kidney tissue

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 49 kDa

Observed band size: 49 kDa