

## ZNF398 Polyclonal Antibody

(Catalog # A63880)

### Background

Function as a transcriptional activator.

### Description

ZNF398 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

Q8TD17

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Zinc finger protein 398 protein (401-642AA)

### Storage

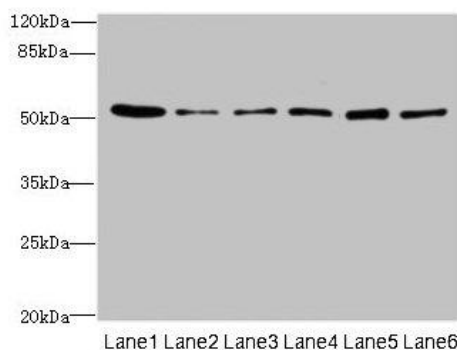
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Zinc finger protein 398, Zinc finger DNA-binding protein p52/p71, ZNF398, KIAA1339, ZER6

### Application

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



#### Western blot

All lanes: ZNF398 antibody at 3ug/ml

Lane 1: 293T whole cell lysate

Lane 2: Hela whole cell lysate

Lane 3: HepG2 whole cell lysate

Lane 4: Jurkat whole cell lysate

Lane 5: Caco-2 whole cell lysate

Lane 6: A549 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 72, 53 kDa

Observed band size: 53 kDa