

---

## ZNF592 Polyclonal Antibody

(Catalog # A52006)

---

### Background

May be involved in transcriptional regulation.

### Description

ZNF592 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

Q92610

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Zinc finger protein 592 protein (1-242AA)

### Storage

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

### Application

ELISA

---