

## polA Polyclonal Antibody

(Catalog # A52038)

### Description

polA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Thermus aquaticus

### Isotype

IgG

### Uniprot ID

P19821

### Purification

Protein G purified

### Immunogen

Recombinant Thermus aquaticus DNA polymerase I, thermostable protein (310-832AA)

### Storage

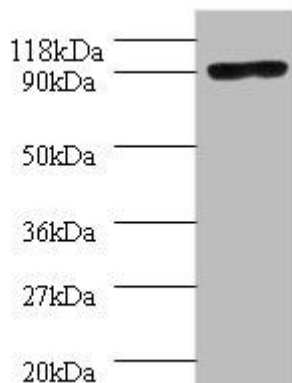
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Taq polymerase 1, polA, pol1

### Application

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



### Western blot

All lanes: polA Polyclonal Antibody at 2ug/ml+293T whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 94kDa

Observed band size: 94kDa