

## APX2 Polyclonal Antibody

(Catalog # A52937)

### Background

Plays a key role in hydrogen peroxide removal.

### Description

APX2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Arabidopsis thaliana

### Isotype

IgG

### Uniprot ID

Q1PER6

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Arabidopsis thaliana L-ascorbate peroxidase 2, cytosolic protein (4-250AA)

### Storage

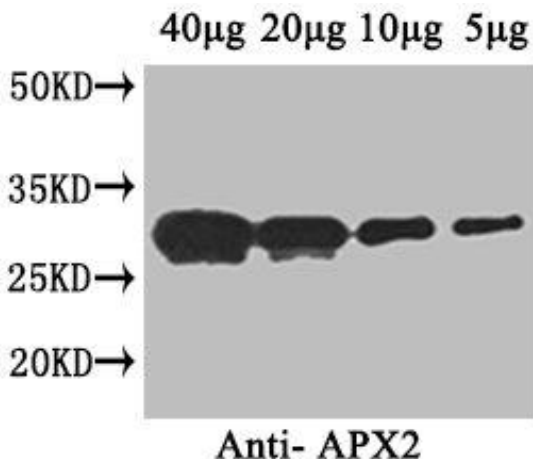
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

L-ascorbate peroxidase 1b, APX2, APX1B, At3g09640, F11F8\_23

### Application

ELISA, WB; Recommended dilution: WB:1:2000-1:5000



### Western Blot

Positive WB detected in: Arabidopsis thaliana (40µg, 20µg, 10µg, 5µg)

All lanes: APX2 Polyclonal Antibody at 1ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 28 kDa

Observed band size: 28 kDa