

## ABI5 Polyclonal Antibody

(Catalog # A54576)

---

### Background

Participates in ABA-regulated gene expression during seed development and subsequent vegetative stage by acting as the major mediator of ABA repression of growth. Binds to the embryo specification element and the ABA-responsive element (ABRE) of the Dc3 gene promoter and to the ABRE of the Em1 and Em6 genes promoters. Can also trans-activate its own promoter, suggesting that it is autoregulated. Plays a role in sugar-mediated senescence.

### Description

ABI5 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

Q9SJN0

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Arabidopsis thaliana Protein ABSCISIC ACID-INSENSITIVE 5 protein (1-442AA)

### 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

Dc3 promoter-binding factor 1 Protein GROWTH-INSENSITIVITY TO ABA 1 bZIP transcription factor 39 ABI5 BZIP39, DPBF1, GIA1, NEM1 At2g36270 F2H17.12

### Application

ELISA

---