

WRKY40 Polyclonal Antibody

(Catalog # A54352)

Background

Transcription factor. Interacts specifically with the W box (5'-(T)TGAC[CT]-3'), a frequently occurring elicitor-responsive cis-acting element

Description

WRKY40 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

Q9SAH7

Purification

>95%, Protein G purified

Immunogen

Recombinant Arabidopsis thaliana Probable WRKY transcription factor 40 protein (1-302AA)

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

WRKY DNA-binding protein 40 WRKY40 At1g80840 F23A5.19

Application

ELISA
