Proteasome 19S Rpn12/S14 subunit polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

Order Online »

BML-PW0440-0025

25µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 26S proteasome non-ATPase regulatory subunit 8

homolog A, AtRPN12a, 26S proteasome regulatory

subunit S14 homolog A

Application WB

Formulation Liquid. Serum.

Host Rabbit

Immunogen Full-length Arabidopsis proteasome 26S (19S regulator

complex) (Rpn12a/S14 subunit).

Species Reactivity Arabidopsis thaliana

Specificity Recognizes the Rpn12a/S14 subunit of the 19S regulator

complex.

UniProt ID Q9SGW3

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

