Proteasome 19S Rpt2/S4 subunit polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

BML-PW0445-0025

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 regulatory subunit 4 homolog A

Application WB

Formulation Liquid. Partially purified antiserum preparation without

sodium azide.

Host Rabbit

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

complex) (Rpt2a/S4 subunit).

Species Reactivity Arabidopsis thaliana

Specificity Recognizes the Rpt2a/S4 subunit of the 19S regulator

complex.

UniProt ID Q9SZD4

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

.

Last modified: May 29, 2024

