# Proteasome activator 11S β subunit polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

# **Ordering Information**

Order Online »

BML-PW8240-0025	25µI
BML-PW8240-0100	100μΙ

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 Proteasome activator complex subunit 2, PA28β

Application IHC, WB

**Formulation** Liquid. Antiserum containing 10mM sodium azide.

GenBank ID 2351198 (mouse)

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 15-31 of mouse

PA28β.

Species Reactivity Human, Mouse, Rat

**Specificity** Recognizes the β subunit of proteasome activator 11S.

UniProt ID P97372 (mouse), Q9UL46 (human)

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

.

Last modified: May 29, 2024

