# Proteasome 19S Rpn2/S1 subunit (human) monoclonal antibody (112-1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

# **Ordering Information**

Order Online »

BML-PW9270-0025	25μΙ
BML-PW9270-0100	100μΙ

Manuals, SDS & CofA

View Online »

# **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name 26S proteasome non-ATPase regulatory subunit 1, 26S

proteasome subunit p112

**Application** WB

**Clone** 112-1

**Formulation** Liquid. In PBS containing 10mM sodium azide.

Gene/Protein Identifier PSMD1 (gene name)

**Host** Mouse

**Immunogen** Recombinant full-length human Rpn2 protein.

lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

**Specificity** Recognizes the Rpn2/S1 subunit of the 19S regulator

complex.

UniProt ID Q99460

Last modified: May 29, 2024



info-

eu@enzolifesciences.com