# Proteasome 20S β1i subunit polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 5

View Online »

# **Ordering Information**

Order Online »

BML-PW8205-0025	25µl
BML-PW8205-0100	100µl

Manuals, SDS & CofA

View Online »

## **Handling & Storage**

Handling Avoid freeze-thaw cycles. Aliquot undiluted antibody into smaller volumes (not less than

 $10\mu L$ ) prior to freezing if appropriate. The use of high quality 'antiserum-grade' plastic or glass vials is recommended. Store diluted antibody at 2-4°C (do not freeze) and use

within 1 month.

Long Term Storage -20°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

**UniProt ID** 

Alternative Name Proteasome subunit β type-9, RING12 protein, LMP2

**Application** IHC, WB

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

Gene/Protein Identifier PSMB9 (gene name)

**Host** Rabbit

Immunogen Recombinant full length mouse proteasome 20S (β1i

subunit).

Species Reactivity Human, Mouse, Rat

**Specificity** Recognizes the β1i subunit of proteasome 20S.

**Technical Info / Product Notes**Various systems for the nomenclature of the proteasome

subunits have been established. This may be a source of confusion as the system on UniProt differs from "standard" nomenclature as described in the literature. The UniProt ID and Gene Name will help to clearly identify the proteins.

P28076 (mouse), P28065 (human)

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com