# Proteasome 20S α3 subunit monoclonal antibody (MCP257)

This antibody is covered by our Worry-Free Guarantee.

Citations: 9

View Online »

# **Ordering Information**

Order Online »

BML-PW8115-0025	25µl
BML-PW8115-0100	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** Proteasome subunit α type-4, Macropain subunit C9,

Proteasome subunit L

**Application** IHC, WB

Clone MCP257

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

**Gene/Protein Identifier** PSMA4 (gene name)

Host Mouse

**Immunogen** Human placenta derived proteasomes.

lgG2b Isotype

**Purity Detail** Partially purified ascites.

Source Purified from hybridoma tissue culture supernatant.

**Species Reactivity** Human, Mouse, Rabbit, Rat

**Specificity** Recognizes the α3 subunit of the 20S proteasome.

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.

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-uk@enzolifesciences.com