

# Proteasome 20S $\beta$ 7 subunit (human) monoclonal antibody (MCP444)

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

BML-PW9150-0100

100 $\mu$ l

## 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  $\beta$  type-4, PROS-26, Macropain  $\beta$  chain

**Application** IP, WB

**Clone** MCP444

**Formulation** Liquid. Ascites in PBS, containing 0.01 M sodium azide.

**Gene/Protein Identifier** PSMB4 (gene name)

**Host** Mouse

**Immunogen** Human placenta derived proteasomes.

**Isotype** IgG1

**Purity Detail** Purified.

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

**Species Reactivity** Human

**Specificity** Recognizes the  $\beta 7$  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.

UniProt ID

P28070

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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