

# Proteasome 20S $\beta$ 2 subunit monoclonal antibody (MCP168)

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.

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

Citations: 15

[View Online »](#)

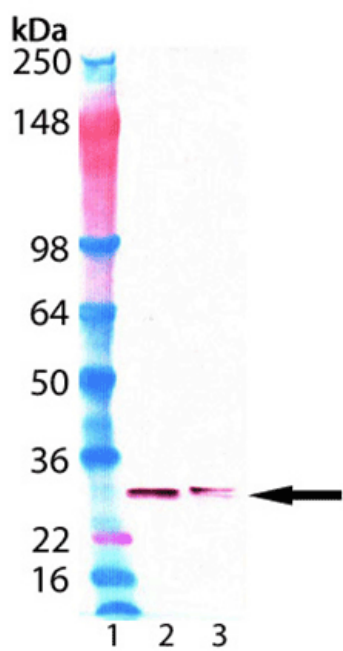
## Ordering Information

[Order Online »](#)

BML-PW8145-0025	25 $\mu$ l
BML-PW8145-0100	100 $\mu$ l

## Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of BML-PW8145: Lane 1: SeeBlue<sup>TM</sup>Plus2 MW Marker, Lane 2: HeLa cell lysate (Prod. No. ADI-LYC-HL100), Lane 3: Jurkat cell lysate (Prod. No. ADI-LYC-JK100).

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Proteasome subunit $\beta$ type-7, Macropain chain Z
Application	IHC, WB
Clone	MCP168
Formulation	Liquid. In PBS containing 10mM sodium azide.
Gene/Protein Identifier	PSMB7 (gene name)
Host	Mouse
Immunogen	Dinitrophenylated proteasomes.
Isotype	IgG1
Purity Detail	Protein G-affinity purified ascites.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Yeast
Specificity	Recognizes the $\beta 2$ subunit of proteasome 20S.
UniProt ID	Q99436

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



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)