

# Proteasome 20S $\alpha$ 2 subunit monoclonal antibody (MCP21)

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

Citations: 6

[View Online »](#)

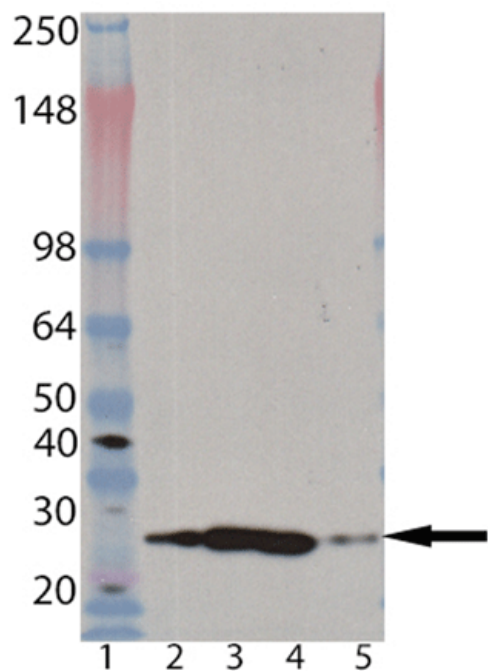
## Ordering Information

[Order Online »](#)

BML-PW8105-0025	25 $\mu$ l
BML-PW8105-0100	100 $\mu$ l

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Proteasome 20S  $\alpha$ 2 subunit, mAb (MCP21) (Prod. No. BML-PW8105): Lane 1: MW marker, Lane 2: HeLa (heat shocked) (Prod. No. ADI-LYC-HL101), Lane 3: Jurkat (Prod. No. ADI-LYC-JK100), Lane 4: HeLa S100 fraction (Prod. No. BML-SW8750), Lane 5: Proteasome 26S (human), (purified) (Prod. No. BML-PW9310).

# 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 $\alpha$ type-2, Macropain subunit C3
Application	IHC, IP, WB
Clone	MCP21
Formulation	Liquid. In PBS containing 10mM sodium azide.
Gene/Protein Identifier	PSMA2 (gene name)
Host	Mouse
Immunogen	Human placenta derived proteasomes.
Isotype	IgG1
Purity Detail	Partially purified ascites.
Source	Purified from mouse ascites.
Species Reactivity	Bovine, Human, Rabbit
Specificity	Recognizes the $\alpha 2$ 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	P25787

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)