

Proteasome 20S β 1 subunit monoclonal antibody (MCP421)

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

Citations: 30

[View Online »](#)

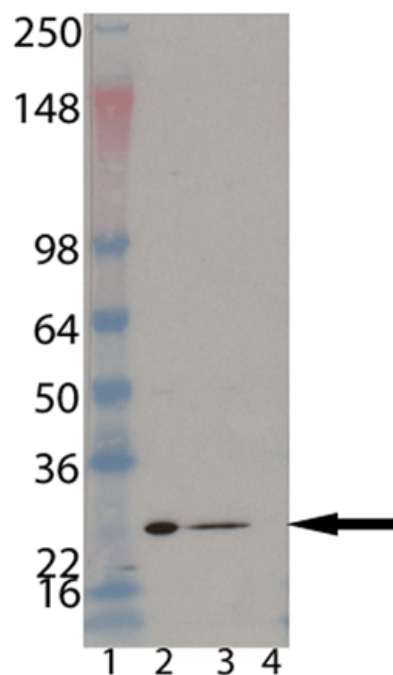
Ordering Information

[Order Online »](#)

BML-PW8140-0025	25 μ l
BML-PW8140-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Proteasome 20S β 1 subunit, mAb (MCP421) (Prod. No. BML-PW8140): Lane 1: MW marker, Lane 2: HeLa (Prod. No. ADI-LYC-HL100), Lane 3: HeLa S100 fraction (Prod. No. BML-SW8750), Lane 4: 3T3 (Prod. No. ADI-LYC-3T100).

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 β type-6, Macropain δ chain, Proteasome subunit Y
Application	IHC, WB
Clone	MCP421
Formulation	Liquid. In PBS containing 10mM sodium azide.
Gene/Protein Identifier	PSMB6 (gene name)
Host	Mouse
Immunogen	Dinitrophenylated proteasomes.
Isotype	IgG1
Purity Detail	Protein G-affinity purified ascites.
Source	Purified from mouse ascites.
Species Reactivity	Human, Rabbit, Yeast
Specificity	Recognizes the β 1 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

P28072

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)