

Proteasome 20S α 6 subunit monoclonal antibody (MCP20)

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

Citations: 20

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW8100-0025	25 μ l
BML-PW8100-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Aliquot undiluted antibody into smaller volumes (not less than 10µL) prior to freezing if appropriate. The use of high quality 'antiserum-grade' plastic or glass vials is recommended. Store diluted antibody at 2-4°C (do not freeze) and use within 1 month. Dilute to working strength with phosphate buffered saline pH 7.2-7.4 and 1% normal goat serum (if a goat anti-mouse IgG linker antibody is to be used).
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PROS30, Proteasome subunit α type-1, Macropain subunit C2, Proteasome γ chain
Application	IHC, IP, WB
Application Notes	<p>Immunoprecipitation: This antibody has been immobilised on CNBr-activated Sepharose CL-4B and used to precipitate proteasomes from HeLa cells. Proteasomal subunits were subsequently separated by 2D PAGE.</p> <p>Western Blot: Single dimension SDS-PAGE of a human placental proteasome preparation or HeLa cell lysate followed by Western blotting gives a single band with a relative molecular weight of approximately 33kDa.</p>
Clone	MCP20
Crossreactivity	Does not cross-react with rat or mouse protein.
Formulation	Liquid. In PBS containing 10mM sodium azide.
Gene/Protein Identifier	PSMA1 (gene name)
Host	Mouse
Immunogen	Human placenta derived proteasomes.
Isotype	IgG1
Purity Detail	Protein G affinity purified.

Source	Purified from ascites.
Species Reactivity	Human, Rabbit
Specificity	Recognizes the $\alpha 6$ subunit of proteasome 20S.
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	P25786
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)