

Proteasome 20S β 2 subunit monoclonal antibody (MCP165)

The hybridoma secreting the antibody to subunit β 2 was generated by fusion of splenocytes from Balb/c mice which had received repeated immunization with dinitrophenylated proteasomes. The antibody (clone MCP165) has been characterised by one- and two dimensional Western blotting. The antibody has not been demonstrated to immunoprecipitate the native 20S proteasome. The immunoglobulin subclass is IgG₁.

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

Citations: 11

[View Online »](#)

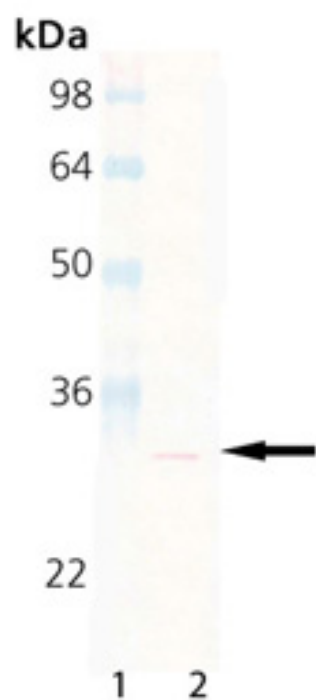
Ordering Information

[Order Online »](#)

BML-PW9300-0025	25 μ l
BML-PW9300-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of Proteasome 20S β 2 subunit, mAb (MCP165), (Prod. No. BML-PW9300): Lane 1: MW Marker, Lane 2: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100)

Handling & Storage

Use/Stability	Store unopened vial at -20°C until required for use. 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.
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 β type-7, Macropain chain Z
Application	IHC, WB
Application Notes	Immunoprecipitation: Does not immunoprecipitate intact proteasomes from HeLa cell preparations. Immunohistochemistry: Has been shown to have applicability in immunohistochemical applications using human tissue. Western blot: Single dimension SDS-PAGE of a HeLa cell lysate followed by Western blotting gives a single band with a relative molecular weight of ~28kDa.
Clone	MCP165
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Gene/Protein Identifier	PSMB7 (gene name)
Host	Mouse
Immunogen	Dinitrophenylated proteasomes.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Mouse
Specificity	Recognizes the β 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	Q99436 (human), P70195 (mouse)
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)