

Proteasome 20S α 4 subunit (human) monoclonal antibody (MCP34)

Two forms of the protein are produced by alternative splicing of the gene - subunits XAPC7-S and XAPC7-L.

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

Citations: 12

[View Online »](#)

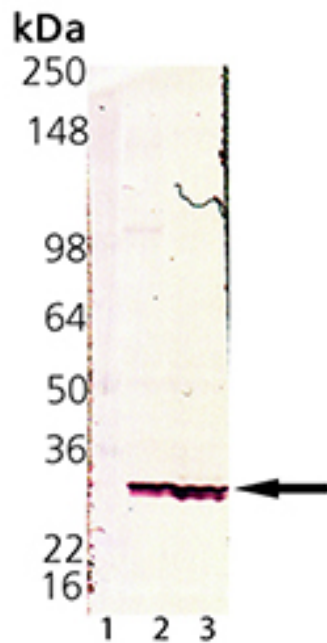
Ordering Information

[Order Online »](#)

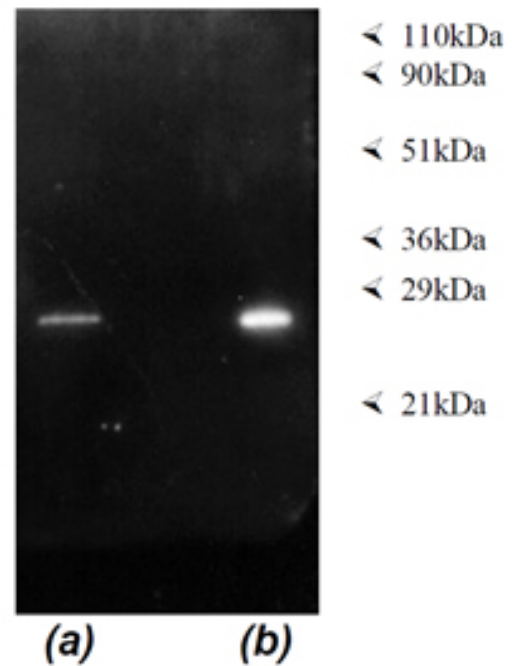
BML-PW8120-0025	25 μ l
BML-PW8120-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of 20S Proteasome subunit α 4, mAb (MCP34) (Prod. No. BML-PW8120): Lane 1: MW marker, Lane 2: HeLa S100 fraction (Prod. No. BML-SW8750), Lane 3: HeLa cell lysate (Prod. No. ADI-LYC-HL100).



Immunoprecipitation analysis: Luminograph of material immunoprecipitated from a HeLa S3 S100 fraction (BML-SW8750) using immobilised BML-PW8335 (lane a) or BML-PW9005 (lane b) after SDS-PAGE followed by blotting onto PVDF and probing with antibody BML-PW8155 to proteasome α and β subunits. Antibody dilution 1:1000 using ECL procedure (1 min exposure).

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).
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Proteasome subunit α type-7, Proteasome subunit RC6-1, Proteasome subunit XAPC7
Application	IHC, IP, WB
Application Notes	<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 28kDa. Upon 2D analysis antibody PW8120 also reacts with two other proteins with the same molecular weight but differing pI. The amount of protein in the spots decreases with their pI. Peptide mapping suggests that the proteins whilst different are closely related. Proteasome preparations containing degradation products often give reaction patterns with several more spots than that shown.</p> <p>Immunoprecipitation: This antibody has been used for immunoprecipitation using standard procedures using immobilised protein A/G for capture. This antibody is also available immobilized on activated agarose (Prod. No. BML-PW9005) and has been used to precipitate proteasomes from a HeLa cell lysate (Prod. No. BML-SW8750).</p>
Clone	MCP34
Formulation	Liquid. In PBS containing 10mM sodium azide.
Gene/Protein Identifier	PSMA7 (gene name)
Host	Mouse
Immunogen	Human placenta derived proteasomes.
Isotype	IgG1

Purity Detail	Partially purified ascites.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from mouse ascites.
Species Reactivity	Human
Specificity	Recognizes the $\alpha 4$ subunit of the 20S proteasome.
Technical Info / Product Notes	<p>The hybridoma secreting the antibody to subunit $\alpha 4$ was generated by fusion of splenocytes from Balb/c mice which had received repeated immunisation with human placenta derived proteasomes. The antibody (clone MCP34) has been extensively characterised by one- and two-dimensional Western blotting and has been demonstrated to immunoprecipitate the native 20S proteasome.</p> <p>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.</p>
UniProt ID	O14818
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)