

Proteasome 20S α 1, 2, 3, 5, 6 & 7 subunits monoclonal antibody (MCP231)

The proteasome is widely recognised as the central enzyme of non-lysosomal protein degradation. It is responsible for intracellular protein turnover and it is also critically involved in many regulatory processes and, in higher eukaryotes, in antigen processing. The 26S proteasome is the key enzyme of the ubiquitin/ATP-dependent pathway of protein degradation. The catalytic core of this unusually large (2000kDa, 450Å in length) complex is formed by the 20S proteasome, a barrel shaped structure shown by electron microscopy to comprise of four rings each containing seven subunits. Based on sequence similarity, all fourteen 20S proteasomal subunit sequences may be classified into two groups, α and β , each group having distinct structural and functional roles. The α -subunits comprise the outer rings and the β -subunits the inner rings of the 20S proteasome. Observations of the eukaryotic proteasome and analysis of subunit sequences indicate that each ring contains seven different subunits (α 7 β 7 β 7 α 7) with a member of each sub-family represented in each particle. Each subunit is located in a unique position within the α - or β -rings. 120S Proteasomes degrade only unfolded proteins in an energy-independent manner, whereas 26S proteasomes degrade native and ubiquitinated proteins in an ATP-dependent manner. The native protein substrates are recognised by subunits, some with ATP binding sites, of the outer 19S caps of the 26S proteasome. The hybridoma secreting the antibody to subunits HC2, HC3, HC8, HC9, Iota and Zeta was generated by fusion of splenocytes from Balb/c mice which had received repeated immunisation with dinitrophenylated proteasomes.

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

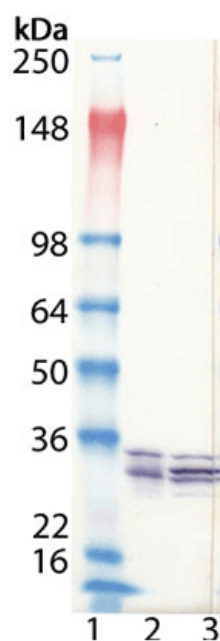
Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS667-0100 | 100µl |
|-----------------|-------|

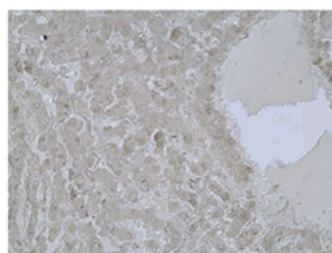
Manuals, SDS & CofA

[View Online »](#)

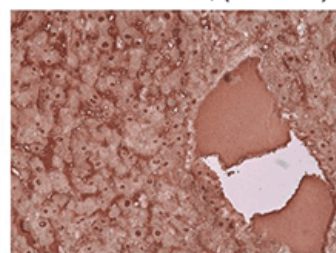


Western Blot analysis of ENZ-ABS667: Lane 1: NovexTMSeeBlueTMPlus2 MW Marker, Lane 2: HeLa cell lysate (Prod. No. ADI-LYC-HL100), Lane 3: Jurkat cell lysate (Prod. No. ADI-LYC-JK100).

Negative Control



Proteasome 20S α 1, 2, 3, 5, 6&7 subunit mAb, (MCP231)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human liver tissue stained with Proteasome 20S α 1, 2, 3, 5, 6 & 7 subunits monoclonal antibody (MCP231), (Prod. No. ENZ-ABS667) at 10 μ g/ml.

Handling & Storage

Handling Store unopened vial at -20°C. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome subunit α type-1, -2, -3, -4, -5 & -6, Macropain subunits C2, C3, C8, C9, ι & ζ

Application IHC, WB

Clone MCP231

Formulation Liquid. In PBS containing 0.01% sodium azide.

Host Mouse

Immunogen Dinitrophenylated proteasomes.

Isotype IgG1 κ

Purity Detail Protein G affinity purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Potato, Rabbit, Rat, Yeast

Specificity Recognizes the α 1, 2, 3, 5, 6 & 7 subunits 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 P25786 (human PSMA1), P25787 (human PSMA2), P25788 (human PSMA3), P25789 (human PSMA4), P28066 (human PSMA5), P60900 (human PSMA6)

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)