

# Proteasome 20S $\beta$ 5i subunit monoclonal antibody (LMP7-1)

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,  $\alpha$  and  $\beta$ , each group having distinct structural and functional roles. The  $\alpha$ -subunits comprise the outer rings and the  $\beta$ -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 ( $\alpha$ 7 $\beta$ 7 $\beta$ 7 $\alpha$ 7) with a member of each sub-family represented in each particle. Each subunit is located in a unique position within the  $\alpha$ - or  $\beta$ -rings. Lmp2, Lmp7 and MECL are interferon gamma-inducible catalytic subunits of the 20S proteasome which may replace the constitutive catalytic subunits, delta, X and Z respectively, during proteasome biogenesis. Lmp2 and Lmp7 alter the cleavage site specificity of the 20S proteasome and are required for the efficient generation of T cell epitopes from a number of viral proteins and for optimal MHC class I cell surface expression .

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

Citations: 21

[View Online »](#)

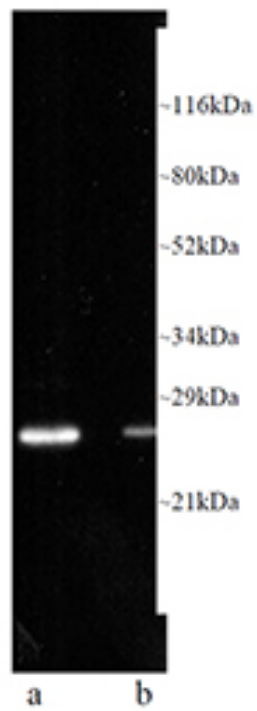
## Ordering Information

[Order Online »](#)

BML-PW8845-0025	25 $\mu$ l
BML-PW8845-0100	100 $\mu$ l

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Luminograph of HeLa cell lysates with (a) and without (b) prior stimulation with  $\gamma$ -interferon after SDS-PAGE followed by blotting onto PVDF and probing with antibody Prod. No. BML-PW8845.

## Handling & Storage

**Long Term Storage**      -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Proteasome subunit $\beta$ type-8, LMP7, RING10 protein, Macropain subunit C13
-------------------------	---

<b>Application</b>	IHC, WB
--------------------	---------

<b>Application Notes</b>	<p>The serum was characterised by single dimension SDS-PAGE using a HeLa cell lysate both with and without prior stimulation with <math>\gamma</math>-interferon. The mature LMP7 protein giving a band at a relative molecular weight of ~25kDa was detected by Western blotting in both stimulated and nonstimulated cells.</p> <p>Not suitable for immunoprecipitation.</p>
--------------------------	--

**Clone** LMP7-1

<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 10mM sodium azide.
--------------------	---

**Gene/Protein Identifier** PSMB8 (gene name), 5696 (Entrez Gene ID)

Host Mouse

<b>Immunogen</b>	Recombinant human proteasome subunit $\beta 5i$ .
------------------	---

**Isotype** IgG1

<b>Purity Detail</b>	Protein G affinity purified.
----------------------	------------------------------

**Source** Purified from ascites.

<b>Species Reactivity</b>	Human, Rat
---------------------------	------------

<b>Specificity</b>	Recognizes the $\beta 5i$ 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

P28062

### 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)