# Proteasome 20S β1i subunit monoclonal antibody (LMP2 – 13)

The proteasome is widely recognised as the central enzyme of nonlysosomal 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 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  $(\alpha7\beta7\beta7\alpha7)$  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: 12

View Online »

### **Ordering Information**

**Order Online** »

BML-PW8840-0025	25µl
BML-PW8840-0100	100µl

Manuals, SDS & CofA

View Online »

# **Handling & Storage**

Long Term Storage -20°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Proteasome subunit β type-9, RING12 protein, LMP2

Application IHC, WB

**Application Notes**The serum was characterised by single dimension SDS-

PAGE using a HeLa cell lysate both with and without prior stimulation with γ-interferon. The mature LMP2 protein of ~22kDa was detected by Western blotting in the γ-

~22kDa was detected by vvestern blotting in the γinterferon stimulated cells and additionally the precursor

form at ~25kDa.

Not suitable for immunoprecipitation.

Clone LMP2-13

**Formulation** Liquid. In PBS containing 10mM sodium azide.

Gene/Protein Identifier PSMB9 (gene name)

**Host** Mouse

**Immunogen** Recombinant human β1i protein.

lgG1

Purity Detail Protein G-affinity purified ascites.

**Source** Purified from ascites.

Species Reactivity Human, Rat

Specificity Recognizes the β1i 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** 

P28065 (human), P28077 (rat)

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-