

Proteasome 19S

Rpt4/S10b subunit

monoclonal antibody

(p42-23)

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 ($\alpha 7\beta 7\beta 7\alpha 7$) with a member of each sub-family represented in each particle. Each subunit is located in a unique position within the α - or β -rings. In addition to the 20S particle, the 26S complex contains over twenty additional proteins, ranging in molecular weight from 25 to 10kDa, located in a distinct complex called the 'PA700 proteasome activator' or the '19S complex', and which determines substrate specificity and provides the multiple enzymatic functions necessary for proteolysis and viability. Systematic analysis of the sub-unit components have revealed at least six members to be ATPases belonging to a new family of ATPbinding proteins, together with a further fifteen sub-units that lack the capacity to bind ATP, isopeptidases and several other proteins thought to be responsible for the unfolding of a protein substrate prior to insertion into the proteolytic core of the 20S proteasome.

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

Citations: 11

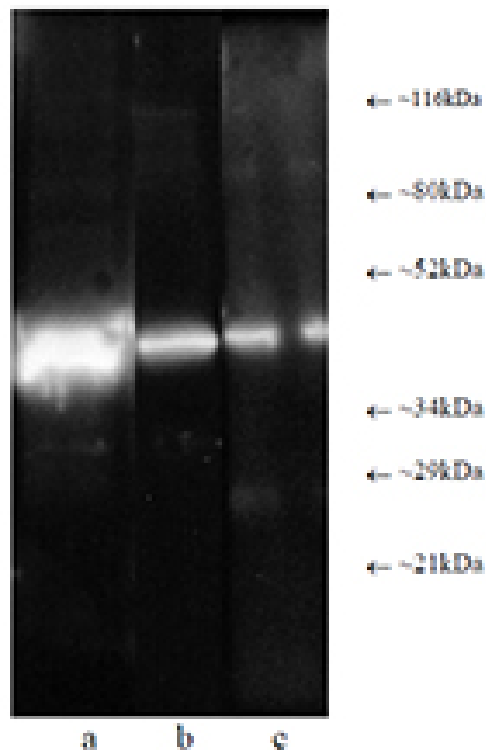
[View Online »](#)

Ordering Information

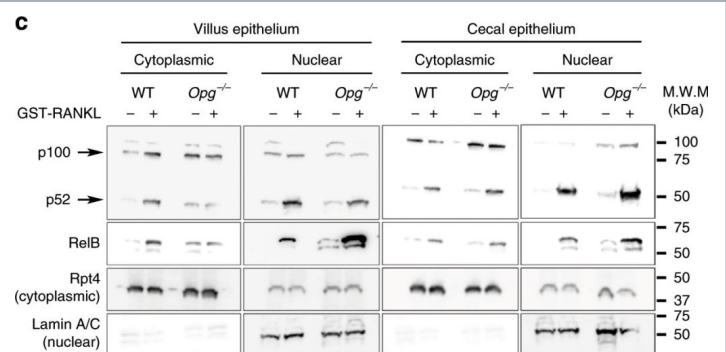
[Order Online »](#)

BML-PW8830-0100

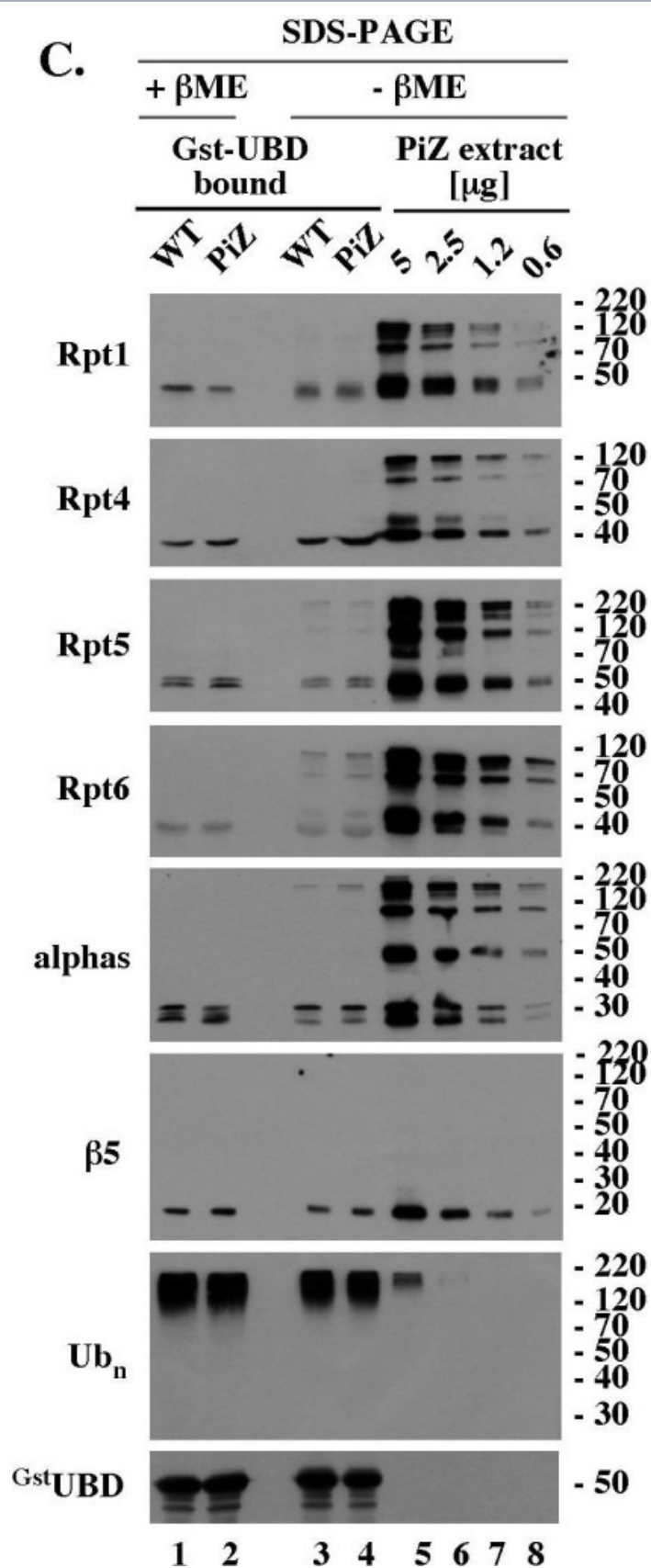
100µl



Western blot analysis: Composite luminograph of (a) HeLa S3 cytosolic preparation, (b) purified 26S proteasome, and (c) human placental proteasome fraction after SDS PAGE followed by blotting onto PVDF membrane and probing with antibody BML-PW8830. Antibody dilution 1:5000 using ECL procedure (1 min exposure).



RANKL–RANK signaling is stimulated in the gut epithelia of *Opg*^{-/-} mice. a GP2⁺ cells (red) are more effectively induced in the cecal epithelium (CE) and ileal villi (VE) of *Opg*^{-/-} mice by RANKL administration. Whole-mount immunohistochemical images of Spi-B (green) and GP2 (red) in the VE and CE of WT (upper panels) and *Opg*^{-/-} mice (lower panels) treated with either GST (control; left) or GST-RANKL (right). Scale bars: 100 μ m. Representative images from three independent experiment are shown. b Quantitative PCR analysis of M-cell marker expression in conventional epithelia from the VE and CE of mice injected with GST (control) or GST-RANKL. Results were normalized to Gapdh expression and are presented relative to the expression in the ileal epithelium from GST-treated mice. Data shown are mean values from three independent experiments (error bars indicate standard deviation). ****p* < 0.005, ***p* < 0.01, **p* < 0.05; *p* values were calculated with the Student's *t*-test (*n* = 3 biologically independent experiments). c, d Nuclear translocation activities of RelB and p52 following RANKL stimulation are enhanced in *Opg*^{-/-} mice. c Western blot analysis of p100/p52 and RelB in the VE and CE of mice injected with GST or GST-RANKL. Rpt4, a subunit of the 26S proteasome, was used as an internal control for the cytoplasmic fraction. Lamin A/C (Lamin) was used as an internal control for the nuclear fraction. Data are representative of two independent experiments. d Right, single confocal planes of the cecal FAE from WT and *Opg*^{-/-} mice. FAE monolayers were stained with anti-RelB (green) and anti-Spi-B (red) antibodies and with Hoechst 33342. Left, a bar graph summarizing the proportions of RelB-positive cells among total numbers of M cells (Spi-B-positive cells). ****p* < 0.005; *p* values were calculated with the Student's *t*-test (*n* = 4 biologically independent experiments). Scale bars: 20 μ m. The source data underlying panels b and d and non-cropped scan images of western blotting (c) are provided as a Source Data file.



Reduction-sensitive modifications typical of aging WT mice accumulate prematurely on selected proteasomal subunits in the livers of PiZ mice. (A). Rpt4 Western blot analysis of unreduced, un-boiled liver extracts. 5 μg of the indicated extracts were mixed with Laemmli buffer without (unreduced samples) or with (reduced samples) β ME (-/+ β ME), separated by SDS-PAGE without prior boiling, and analyzed by Western blot with antibodies specific to Rpt4. (B). Rpt4 Western blot analysis of unreduced, but boiled, samples. Experiment like in A, lanes 1–9, except that extracts were mixed with Laemmli buffer without β ME (- β ME) and boiled for 4

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome 26S subunit ATPase 6, Proteasome subunit p42, 26S protease regulatory subunit 10B

Application WB

Application Notes Single dimension SDS-PAGE of proteasomal preparations from a variety of sources followed by Western blotting gives a single band with a relative molecular weight of approximately 44kDa.
Not suitable for immunoprecipitation.

Clone p42-23

Formulation Liquid. Contains 10mM sodium azide.

Gene/Protein Identifier PSMC6 (gene name)

Host Mouse

Immunogen Recombinant human Rpt4.

Isotype IgG2a

Purity Detail Partially purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse

Specificity Recognizes the Rpt4/S10b subunit of the 19S regulator complex.



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.