

# Proteasome activator

## 11S $\alpha$ subunit

### polyclonal antibody

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. An enzymatic cascade is responsible for the attachment of multiple ubiquitin molecules to lysine residues of proteins targeted for degradation. 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 (31 subunits) is formed by the 20S proteasome, a barrel shaped structure shown by electron microscopy to comprise of four rings each containing seven subunits. 20S Proteasomes degrade only unfolded proteins in an energy-independent manner, whereas 26S proteasomes degrade native and ubiquitylated 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 19S regulatory subunit recognizes ubiquitinated proteins and plays an essential role in unfolding and translocating targets into the lumen of the 20S subunit. A second activator which can associate with the 20S proteasome in the absence of ATP is known as PA28 or the 11S regulator. The pure PA28 activator is a complex of two alternating subunits, PA28 $\alpha$  and PA28 $\beta$ , which share approximately 50% homology but also show considerable similarity (30-40%) to a nuclear protein of unknown function, the Ki autoantigen (recently referred to as PA28 $\gamma$ ). These subunits, with an apparent relative molecular weight of approximately 29kDa, form ringlike heteromeric complexes of ~200kDa possibly with an  $\alpha_3\beta_3$  stoichiometry. Electron microscopic studies have shown PA28 to be a ring shaped particle which, like the 19S, caps the 20S proteasome, by binding to the  $\alpha$ -rings, at both or either end. The complex may, however, be readily dissociated. The finding that PA28 modulates the proteasome-catalysed production of antigenic peptides presented to the immune system on MHC class I molecules indicates a cellular function of this activator in antigen processing. Several genetic diseases are associated with defects in the ubiquitin-proteasome pathway. Some examples of affected proteins include those linked to cystic fibrosis, Angelman's syndrome, and Liddle syndrome.

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

Citations: 15

[View Online »](#)

Ordering Information

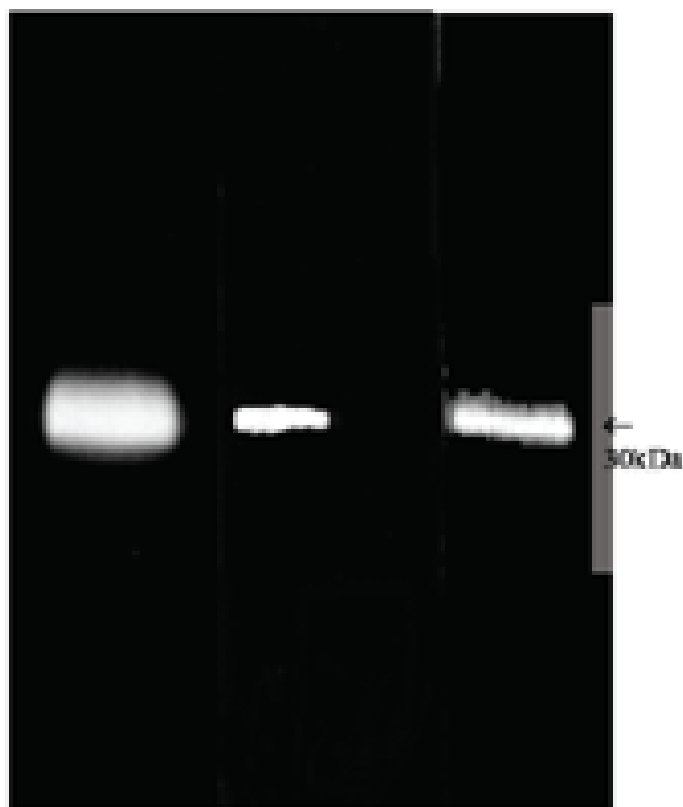
[Order Online »](#)

BML-PW8185-0025	25µl
BML-PW8185-0100	100µl

Manuals, SDS & CofA

[View Online »](#)





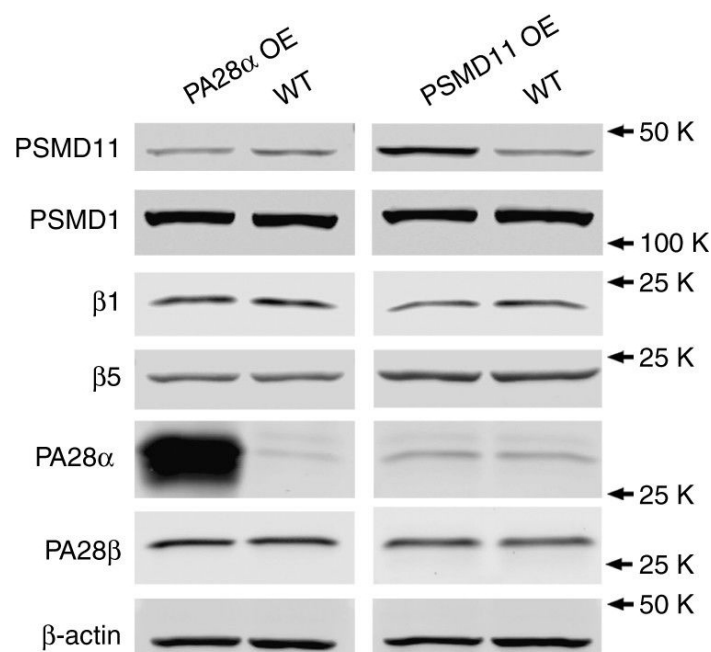
(a)

(b)

(c)

Luminograph of (a) mouse liver preparation, (b) human placental proteasome preparation, and (c) HeLa cell lysate after PAGE followed by blotting onto nitrocellulose and probing with antibody BML-PW8185. Antibody dilution 1:1000 using ECL procedure (1 min exposure).

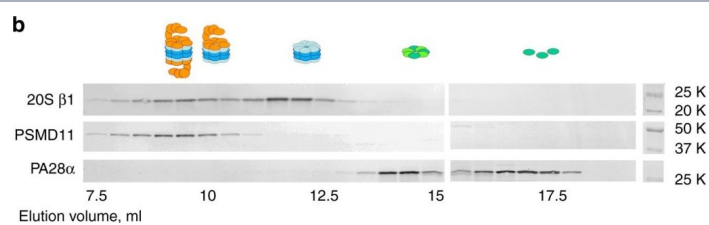
a



#### Characterization of PA28α and PSMD11

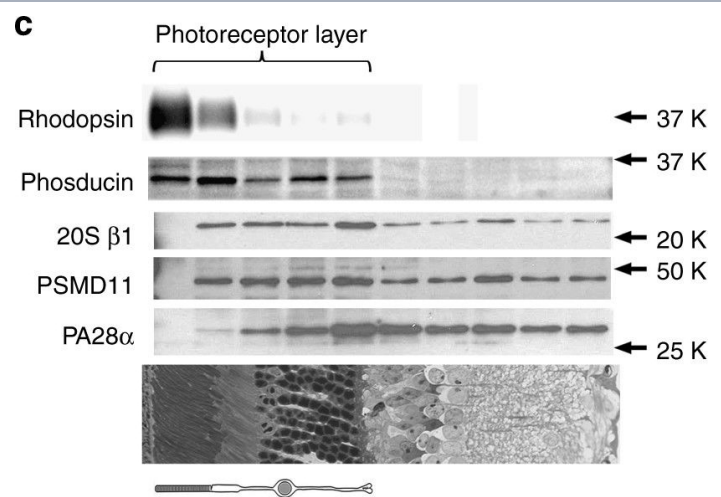
overexpressing (OE) mice. a Western blots of proteasomal subunits in retinal lysates containing 30  $\mu$ g total protein. Bands were visualized using the LiCor Odyssey imaging system. Each protein was analyzed in at least 3 pairs of 1-month-old WT and overexpressing animals. b Retinal morphology of 3-month-old overexpressing and WT mice. Retinas were embedded in plastic, 1  $\mu$ m cross-sections were stained by toluidine blue and analyzed by light microscopy. Data are taken from one of the five similar experiments; scale bar: 20  $\mu$ m. c Chymotrypsin-like proteasomal activity in retinal extracts from 1-month-old overexpressing and WT mice; measurements were performed in the presence or absence of ATP, as indicated. The number of measurements was 10, 7, and 5 for WT, PA28α overexpressing, and PSMD11 overexpressing mice, respectively. The data are shown as mean  $\pm$  SEM; p values determined across individual preparations are indicated in the text. d Fractionation of proteasomal components in retinal extracts from 2-month-old overexpressing and WT mice by size-exclusion chromatography on a Superose-6 Increase column. Proteins in 0.5 ml fractions were probed by western blotting using antibodies against the  $\beta$ 1 subunit of the 20S proteasome core, PSMD11 subunit of the 19S proteasome cap, and PA28α subunit of the 11S cap. Data are taken from one of the three similar experiments

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Increased proteasomal activity supports photoreceptor survival in inherited retinal degeneration. *Nat Commun* (2018)



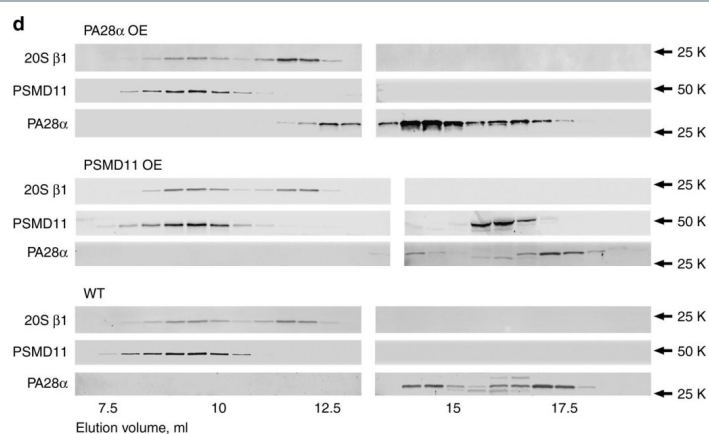
**b** Proteasome composition of the mouse retina. **a** The molar ratio among 20S, 19S, and 11S proteasomal components determined by quantitative mass spectrometry. Data are shown as mean  $\pm$  SEM;  $n = 3$ . **b** Fractionation of proteasome components in retinal extracts from 1-month-old mice (200  $\mu$ g total protein) by size-exclusion chromatography on a Superose-6 column. Proteins in 0.5 ml fractions were probed by western blotting using antibodies against  $\beta 1$  subunit of the 20S proteasome core, PSMD11 subunit of the 19S proteasome cap, and PA28 $\alpha$  subunit of the 11S cap. Data are taken from one of the four similar experiments. **c** The distribution of  $\beta 1$ , PSMD11, and PA28 $\alpha$  in 20  $\mu$ m serial tangential sections throughout the entire WT mouse retina. Each section was solubilized in 30  $\mu$ l SDS-PAGE sample buffer for analysis. Proteins were visualized by western blotting using the ECL technique. Rhodopsin was used as a photoreceptor outer segment marker; phosducin was used as a marker of the entire photoreceptor layer. Data are taken from one of two similar experiments. A representative retinal cross-section is shown below western blot panes; the corresponding position of the photoreceptor cells is illustrated by a cartoon

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Increased proteasomal activity supports photoreceptor survival in inherited retinal degeneration. *Nat Commun* (2018)



**c** Proteasome composition of the mouse retina. **a** The molar ratio among 20S, 19S, and 11S proteasomal components determined by quantitative mass spectrometry. Data are shown as mean  $\pm$  SEM;  $n = 3$ . **b** Fractionation of proteasome components in retinal extracts from 1-month-old mice (200  $\mu$ g total protein) by size-exclusion chromatography on a Superose-6 column. Proteins in 0.5 ml fractions were probed by western blotting using antibodies against  $\beta 1$  subunit of the 20S proteasome core, PSMD11 subunit of the 19S proteasome cap, and PA28 $\alpha$  subunit of the 11S cap. Data are taken from one of the four similar experiments. **c** The distribution of  $\beta 1$ , PSMD11, and PA28 $\alpha$  in 20  $\mu$ m serial tangential sections throughout the entire WT mouse retina. Each section was solubilized in 30  $\mu$ l SDS-PAGE sample buffer for analysis. Proteins were visualized by western blotting using the ECL technique. Rhodopsin was used as a photoreceptor outer segment marker; phosducin was used as a marker of the entire photoreceptor layer. Data are taken from one of two similar experiments. A representative retinal cross-section is shown below western blot panes; the corresponding position of the photoreceptor cells is illustrated by a cartoon

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Increased proteasomal activity supports photoreceptor survival in inherited retinal degeneration. *Nat Commun* (2018)



## Characterization of PA28 $\alpha$ and PSMD11

overexpressing (OE) mice. a Western blots of proteasomal subunits in retinal lysates containing 30  $\mu$ g total protein. Bands were visualized using the LiCor Odyssey imaging system. Each protein was analyzed in at least 3 pairs of 1-month-old WT and overexpressing animals. b Retinal morphology of 3-month-old overexpressing and WT mice. Retinas were embedded in plastic, 1  $\mu$ m cross-sections were stained by toluidine blue and analyzed by light microscopy. Data are taken from one of the five similar experiments; scale bar: 20  $\mu$ m. c Chymotrypsin-like proteasomal activity in retinal extracts from 1-month-old overexpressing and WT mice; measurements were performed in the presence or absence of ATP, as indicated. The number of measurements was 10, 7, and 5 for WT, PA28 $\alpha$  overexpressing, and PSMD11 overexpressing mice, respectively. The data are shown as mean  $\pm$  SEM; p values determined across individual preparations are indicated in the text. d Fractionation of proteasomal components in retinal extracts from 2-month-old overexpressing and WT mice by size-exclusion chromatography on a Superose-6 Increase column. Proteins in 0.5 ml fractions were probed by western blotting using antibodies against the  $\beta$ 1 subunit of the 20S proteasome core, PSMD11 subunit of the 19S proteasome cap, and PA28 $\alpha$  subunit of the 11S cap. Data are taken from one of the three similar experiments

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Increased proteasomal activity supports photoreceptor survival in inherited retinal degeneration. *Nat Commun* (2018)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** PA28α, Proteasome activator complex subunit 1

**Application** WB

**Formulation** Liquid. Antiserum containing 10mM sodium azide.

**GenBank ID** 1698570 (mouse)

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 5-19 of mouse PA28α.

**Species Reactivity** Human, Mouse, Rat

**Specificity** Recognizes the α subunit of proteasome activator 11S.

**UniProt ID** P97371 (mouse), Q06323 (human)

**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-  
fr@enzolifesciences.com](mailto:info-fr@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)