

# Proteasome 19S

## Rpn10/S5a subunit

### monoclonal antibody

### (S5a-18)

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

Citations: 13

[View Online »](#)

#### Ordering Information

[Order Online »](#)

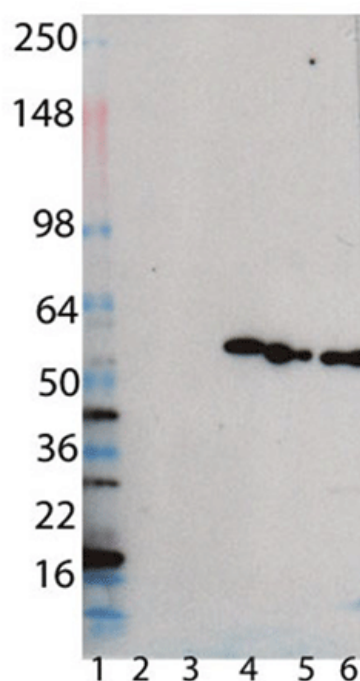
BML-PW9250-0025	25µl
BML-PW9250-0100	100µl

Manuals, SDS & CofA

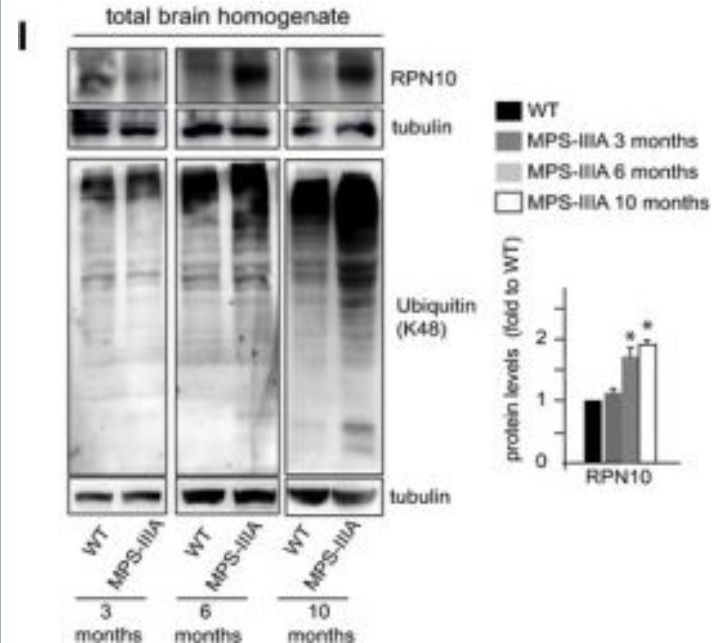
[View Online »](#)







**Figure 1.** Western blot analysis of Proteasome 19S Rpn10/S5a subunit, mAb (S5a-18) (Prod. No. BML-PW9250): Lane 1: MW marker, Lane 2: 3T3 (Prod. No. ADI-LYC-3T100), Lane 3: L929, Lane 4: PC-12 (Prod. No. ADI-LYC-PC100), Lane 5: MDBK, Lane 6: HeLa S100 fraction (Prod. No. BML-SW8750).



Lysosomal driven deregulation of  $\alpha$  synuclein and CSP $\alpha$  degradation WB analysis of LC3 and p62 (an autophagy substrate) was performed on WT and MPS IIIA brain samples at the indicated ages. WB quantitation is shown.  $\alpha$  Synuclein was immunoblotted in WT and MPS IIIA in both total and synaptosomal brain fractions at the indicated ages after sequential extraction with detergents with increased strength. Soluble (Sol.), lowly insoluble (L. Insol.), and highly insoluble (H. Insol.) forms correspond to the protein solubilized in Triton X 100, 10% SDS and 8 M urea, respectively. Co immunofluorescence analysis of  $\alpha$  synuclein with SMI 32 in WT and MPS IIIA hippocampal neurons (DIV14).  $\alpha$  Synuclein synaptic puncta present in a neurite tract of 10  $\mu$ m is shown in a representative enlarged image. Quantification of  $\alpha$  synuclein synaptic puncta was calculated from 30 different enlarged images. Confocal analysis of  $\alpha$  synuclein (green) and LAMP1 (red) in WT and MPS IIIA hippocampal neurons (DIV14). Enlarged merge images are also shown. Co localization was quantified using the MCC (ImageJ) and displayed as percentage ( $MCC \times 100$ ) of  $\alpha$  synuclein co localizing with LAMP1 (15 different images taken from 4 to 5 coverslips for each group). CSP $\alpha$  was immunoblotted in WT and MPS IIIA total brain lysates at the indicated ages after sequential extraction with detergents with increased strength as in (B). Co immunofluorescence analysis of CSP $\alpha$  and SMI 32 in DIV14 hippocampal neurons. CSP $\alpha$  synaptic puncta was quantified as in (C). CSP $\alpha$  protein levels were evaluated by immunoblot analysis in WT and MPS IIIA hippocampal neurons (DIV14) at different times after cycloheximide treatment and expressed as percentage of remaining protein at T0 (100%). The proteasome was inhibited as indicated. WB quantification is shown. Palmitoylation dependent shift in



## Handling & Storage

**Handling** Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 26S proteasome non-ATPase regulatory subunit 4, Antisecretory factor 1, Multiubiquitin chain-binding protein

**Application** IP, WB

**Clone** S5a-18

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

**Gene/Protein Identifier** PSMD4 (gene name)

**Host** Mouse

**Immunogen** Recombinant full-length human Rpn10 protein.

**Isotype** IgG1

**Recommendation Dilutions/Conditions** Western Blot (1:1,000, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from hybridoma tissue culture supernatant.

**Species Reactivity** Bovine, Human, Rat

**Specificity** Recognizes the Rpn10/S5a subunit of the 19S regulator complex.

**UniProt ID** P55036



ENZO LIFE SCIENCES,  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 1 42 44 55 55  
[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)

**This antibody is covered by our Worry Free Guarantee**