

Proteasome 19S Rpt3/S6b subunit monoclonal antibody (TBP7-27)

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

Citations: 3

[View Online »](#)

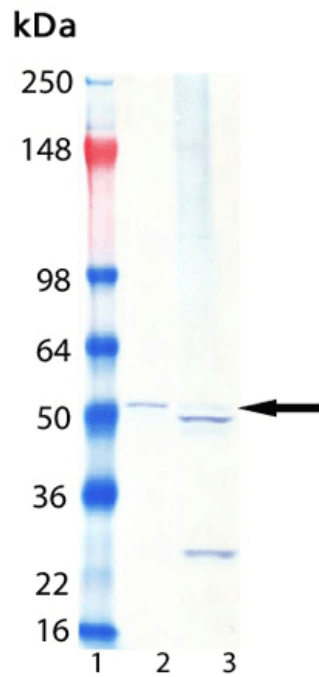
Ordering Information

[Order Online »](#)

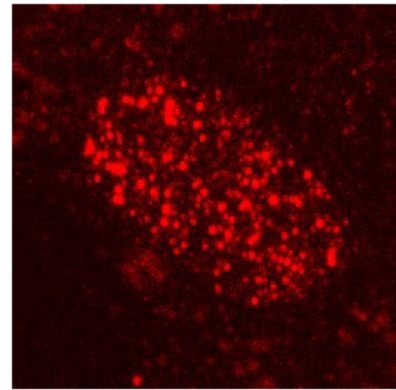
BML-PW8765-0025	25µl
BML-PW8765-0100	100µl

Manuals, SDS & CofA

[View Online »](#)



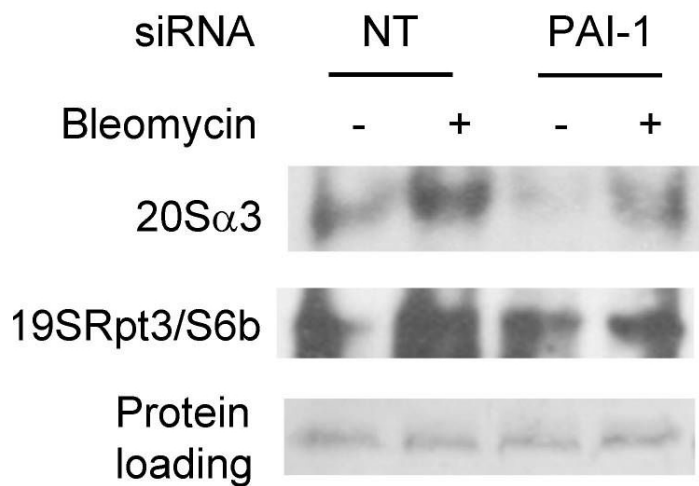
Western Blot analysis of BML-PW8765: Lane 1: MW Marker, Lane 2: HeLa lysate, (Prod. No. ADI-LYC-HL100), Lane 3: HeLa S100 fraction, (Prod. No. BML-SW8750).



This antibody has been used for immunofluorescence. The image was captured on a confocal microscope and is a projection of 17 images taken at 0.4 μm intervals through the cell nucleus. The cells are human primary fibroblasts fixed in 4% paraformaldehyde and permeabilised in 0.5% Triton X-100. BML-PW8765 was used at a dilution of 1:1000 and a goat anti-mouse secondary antibody labeled with fluorophore 594 (Molecular Probes). (Data courtesy of Dr. Carol Shiels, Imperial College, London.)

A

IP with PAI-1 Ab

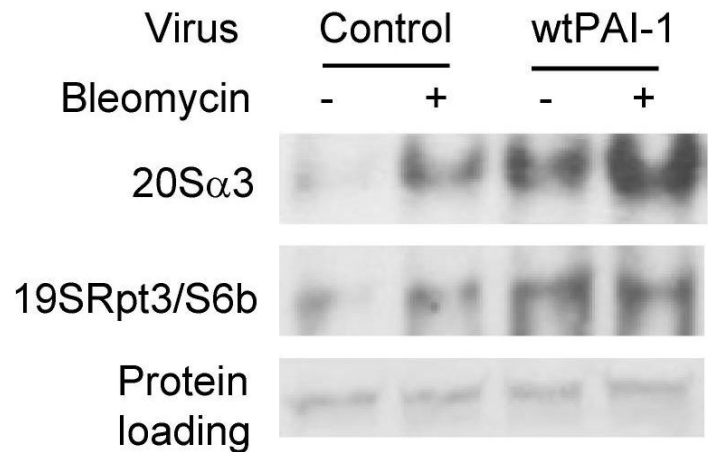


PAI-1 binds to proteasome components in A549 cells. (A,D) Immunoprecipitation-immunoblotting analysis of PAI-1 interaction with proteasome 20S α 3 and 19S Rpt3/S6b subunits in A549 cells. A549 cells were transfected with PAI-1 siRNA/non-target (NT) siRNA or transduced with control/PAI-1 expressing viruses and then treated with bleomycin. PAI-1 protein was immunoprecipitated with anti-mouse PAI-1 monoclonal antibody, and Westerns were conducted with specific antibody to 20S α 3 or 19SRpt3/S6b. Proteins on the membrane were stained with Ponceau S to show equal sample loading. (B,C,E,F) Semi-quantification of the band intensities, normalized by the corresponding protein staining band (n = 3).

Image collected and cropped by CiteAb under a CC-BY license from the following publication: PAI-1 Regulation of p53 Expression and Senescence in Type II Alveolar Epithelial Cells. *Cells* (2023)

D

IP with PAI-1 Ab



PAI-1 binds to proteasome components in A549 cells. (A,D) Immunoprecipitation-immunoblotting analysis of PAI-1 interaction with proteasome 20S α 3 and 19S Rpt3/S6b subunits in A549 cells. A549 cells were transfected with PAI-1 siRNA/non-target (NT) siRNA or transduced with control/PAI-1 expressing viruses and then treated with bleomycin. PAI-1 protein was immunoprecipitated with anti-mouse PAI-1 monoclonal antibody, and Westerns were conducted with specific antibody to 20S α 3 or 19SRpt3/S6b. Proteins on the membrane were stained with Ponceau S to show equal sample loading. (B,C,E,F) Semi-quantification of the band intensities, normalized by the corresponding protein staining band (n = 3).

Image collected and cropped by CiteAb under a CC-BY license from the following publication: PAI-1 Regulation of p53 Expression and Senescence in Type II Alveolar Epithelial Cells. *Cells* (2023)

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 26S protease regulatory subunit 6B, Proteasome 26S subunit ATPase 4, Tat-binding protein 7, TBP7

Application IF, IHC, IP, WB

Clone TBP7-27

Formulation Liquid. Contains 10mM sodium azide.

Host Mouse

Immunogen Recombinant full-length human Rpt3.

Isotype IgG1

Purity Detail Partially purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Rabbit

Specificity Recognizes the Rpt3/S6b subunit of the 19S regulator complex.

UniProt ID P43686

Worry-free Guarantee

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



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-
be@enzolifesciences.com](mailto:info-be@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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.