

Proteasome 19S

Rpt2/S4 subunit

(human) polyclonal

antibody

The 19S regulatory complex consists of a 'base' sub-complex, composed of a hetero-hexameric ring of AAA-ATPases that not only function as the gate-opener to the 20S proteasome but also, in conjunction with two other high-molecular-mass non-ATPase subunits (Rpn2/S1 and Rpn1/S2), as a platform for protein unfolding. The 'lid' sub-complex of the 19S complex, composed of non-ATPase subunits is involved in the binding and deubiquitinylation of proteasome targeted ubiquitin conjugated proteins and assists in the transport of substrates through the 'base' and into the 20S catalytic core.

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

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW0530-0025	25µl
BML-PW0530-0100	100µl

Manuals, SDS & CofA

[View Online »](#)

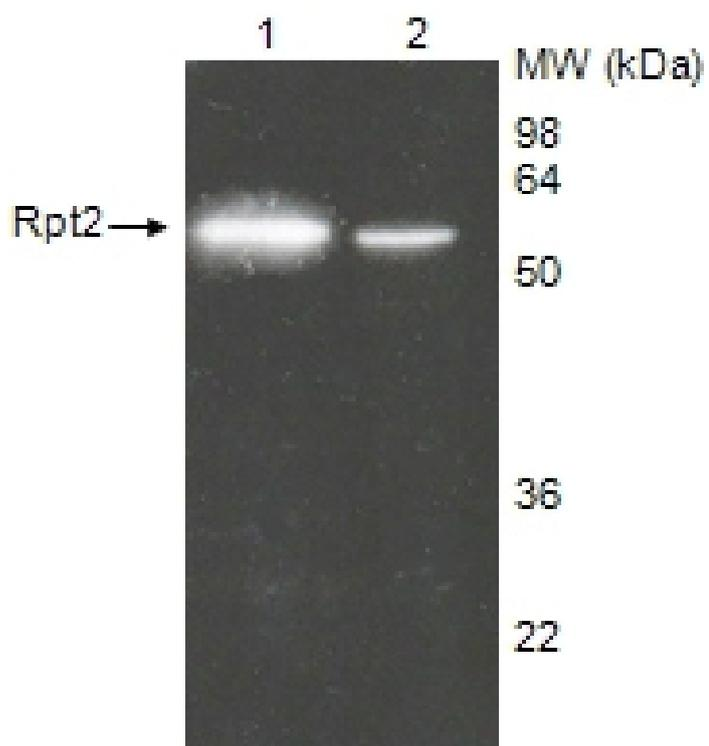


Figure: Western blot showing detection of Rpt2/S4 19S regulatory subunit. **Lane 1:** Purified 26S proteasome (Prod. No. [BML-PW9310](#), 1 μ g). **Lane 2:** HeLa S100 cytosolic fraction (Prod. No. BML-SW8750)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C. Dilute with PBS, pH 7.2-7.4, containing 1% normal goat serum (if a goat anti-rabbit IgG linker antibody is to be used). Store diluted antibody at +4°C, do not freeze and use within 1 month.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Proteasome 26S subunit ATPase 1, 26S protease regulatory subunit 4
Application	WB
Formulation	Liquid. In PBS containing 0.01% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 368-383 of human proteasome 19S (Rpt2/S4 subunit).
Purity Detail	Partially purified by salt precipitation.
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
UniProt ID	P62191
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-
be@enzolifesciences.com](mailto:info-be@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)