

Proteasome 19S Rpn12/S14 subunit (human) monoclonal antibody (P31-27)

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

Manuals, SDS & CofA

[View Online »](#)

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 proteasome non-ATPase regulatory subunit 8

Application IHC, WB

Clone p31-27

Formulation Liquid. Containing 10mM sodium azide.

Gene/Protein Identifier PSMD8 (gene name)

Host Mouse

Immunogen Recombinant Rpn12 protein.

Isotype IgG1

Purity Detail Partially purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

Specificity Recognizes human proteasome 26S (19S regulator complex) (Rpn12/S14 subunit).

UniProt ID P48556

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)