

Proteasome 19S Rpt1/S7 subunit (human) monoclonal antibody (MSS1-92)

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

Citations: 1

[View Online »](#)

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 protease regulatory subunit 7, Proteasome 26S subunit ATPase 2, MSS1

Application IP

Clone MSS1-92

Formulation Liquid. In TRIS buffer containing 10mM sodium azide.

Host Mouse

Immunogen Recombinant human Rpt1 protein.

Isotype IgG1

Purity Detail Ammonium sulphate precipitation and ion-exchanging chromatography purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

Specificity Recognizes the Rpt1/S7 subunit of the 19S regulator complex.

UniProt ID P35998

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)