

Proteasome 20S (alpha subunits) antibody sampler pack

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents

Contains 10 µl each of:

Proteasome 20S α6 subunit, mAb (MCP20) (Prod. No. [BML-PW8100](#))

Proteasome 20S α2 subunit, mAb (MCP21) (Prod. No. [BML-PW8105](#))

Proteasome 20S α7 subunit, mAb (MCP72) (Prod. No. [BML-PW8110](#))

Proteasome 20S α3 subunit, mAb (MCP257) (Prod. No. [BML-PW8115](#))

Proteasome 20S α4 subunit (human), mAb (MCP34) (Prod. No. [BML-PW8120](#))

Proteasome 20S α5 subunit, mAb (MCP196) (Prod. No. [BML-PW8125](#))

Proteasome 20S α1, 2, 3, 5, 6 & 7 subunits, mAb (MCP231) (Prod. No. [BML-PW8195](#))

Proteasome 20S core subunits, pAb (Prod. No. [BML-PW8155](#))

Source

Provided in individual antibody datasheet.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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