

Proteasome 20S β 1i subunit polyclonal antibody

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------------|
| BML-PW8205-0025 | 25 μ l |
| BML-PW8205-0100 | 100 μ l |

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|-------------------|---|
| Handling | Avoid freeze-thaw cycles. Aliquot undiluted antibody into smaller volumes (not less than 10µL) prior to freezing if appropriate. The use of high quality 'antiserum-grade' plastic or glass vials is recommended. Store diluted antibody at 2-4°C (do not freeze) and use within 1 month. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|--------------------------------|---|
| Alternative Name | Proteasome subunit β type-9, RING12 protein, LMP2 |
| Application | IHC, WB |
| Formulation | Liquid. Antiserum containing 10mM sodium azide. |
| Gene/Protein Identifier | PSMB9 (gene name) |
| Host | Rabbit |
| Immunogen | Recombinant full length mouse proteasome 20S (β 1i subunit). |
| Species Reactivity | Human, Mouse, Rat |
| Specificity | Recognizes the β 1i subunit of proteasome 20S. |
| Technical Info / Product Notes | Various systems for the nomenclature of the proteasome subunits have been established. This may be a source of confusion as the system on UniProt differs from "standard" nomenclature as described in the literature. The UniProt ID and Gene Name will help to clearly identify the proteins. |
| UniProt ID | P28076 (mouse), P28065 (human) |

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



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