

POMP polyclonal antibody

The assembly of individual mammalian proteasome subunits into catalytically active 20S proteasome is a complex process, with biogenesis of the mammalian 20S proteasome occurring *via* precursor complexes containing α and unprocessed β subunits. The pro-peptides of the β subunits are thought to prevent premature conversion of the precursor complexes into matured particles and are needed for efficient β subunit incorporation. The complex biogenesis is tightly regulated, requiring additional components such as the maturation factor POMP (proteasome maturation protein), an ubiquitous protein in eukaryotic cells. Mammalian POMP is functionally related to but not interchangeable with its yeast homologue Ump1. POMP is associated purely with the precursor intermediates and is degraded upon final maturation. Northern blot experiments have revealed that expression of POMP is induced after treatment with interferon γ .

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

Ordering Information

[Order Online »](#)

BML-PW9715-0025	25 μ l
BML-PW9715-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome maturation protein, UMP1 homolog

Application IHC, IP, WB

Formulation Liquid. In PBS containing 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 129-141 of human POMP.

Purity Detail Purified by salt precipitation and peptide immunochromatography.

Species Reactivity Human, Monkey, Rat

UniProt ID Q9Y244

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



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)