

Proteasome 20S (human), (purified)

High integrity preparation for use in
proteasome research.

The 20S proteasome is the catalytic core of the 26S proteasome, the central protease of the ubiquitin pathway of protein degradation. Unlike the 26S proteasome, which degrades proteins in an ATP-dependent manner, the 20S proteasome is ATP-independent and only degrades entirely unfolded polypeptides.

Citations: 32

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW8720-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. 20S Proteasome remains relatively stable and active so long as repetitive freeze-thaw cycles are avoided. It has been demonstrated that 20S proteasome undergoes only slow loss of activity when stored at 4°C for up to six months. If it is to be used on a regular basis, it is recommended that the material should be stored between 0°C-4°C during the duration of use. For longer term storage the material may be dispensed as single use aliquots and stored at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Concentration	1.0mg/ml
Formulation	Liquid. In 20mM TRIS-HCl, pH 7.2, containing 1mM EDTA, 1mM DTT, 1mM sodium azide and 50% glycerol.
Source	Isolated and purified from human erythrocytes.



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)