

Proteasome 26S (human), (purified)

High integrity preparation for use in
proteasome research.

Highly purified preparation of '26S' proteasomes useful for carrying out *in vitro* protein degradation studies with suitably ubiquitinated protein substrates.

Citations: 11

[View Online »](#)

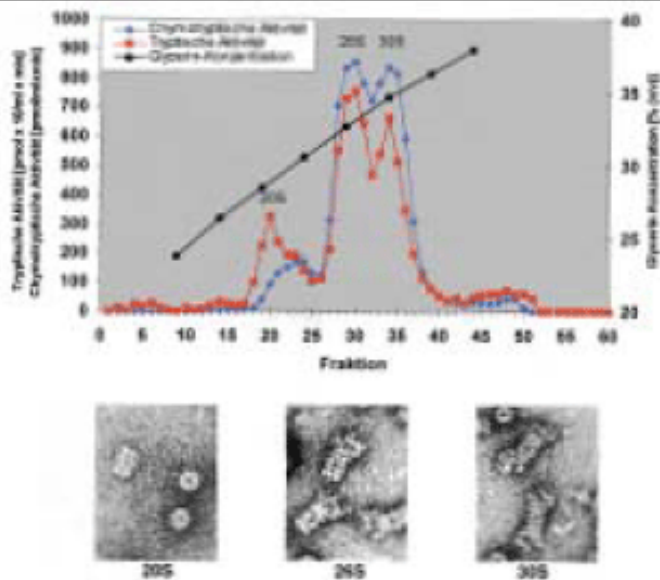
Ordering Information

[Order Online »](#)

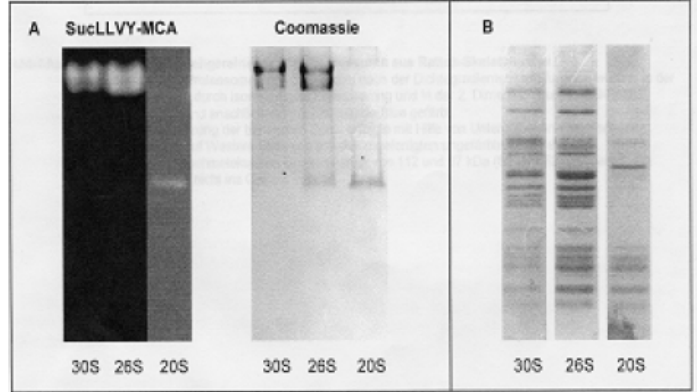
BML-PW9310-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



A typical example of a sucrose density gradient separation and purification of 20S, 26S and 30S proteasome derived from human erythrocytes together with electron micrographs of the individual complexes. The protein concentration used for micrography is best at ~2.5µg/mL.



A – Typical results from substrate overlay (carried out in buffer containing ATP and an ATP-regenerating system) and Coomassie staining of non denaturing gel of products isolated according to the purification shown alongside. Protein concentration is ~15-20µg per lane..
B – Coomassie staining of denaturing gel showing presence of proteasomal subunits.

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. When ready for use the enzyme should be thawed by standing on ice. If the enzymatic activity of the 26S proteasome is to be measured, it should be used immediately after thawing since the enzyme complex is labile. After dissociation of the 26S complex the 20S proteasome activity is relatively stable.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation	Suspended in TSD buffer (50µg in 10mM TRIS, containing 25mM potassium chloride, 1.1mM magnesium chloride, 0.1mM ethylenediaminetetraacetic acid, 1mM dithiothreitol, 1mM sodium azide, 2mM ATP, pH 7.0, and 35% glycerol).
Source	Isolated from human erythrocytes. Consists of a high purity mixture of '26S' proteasomes singly (26S) and doubly (30S) capped with 19S regulatory subunit complexes in the ratio of 40% single cap : 60% double capped at the time of preparation.

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)