

Immunoproteasome

20S (human), (purified)

The eukaryotic proteasome contains 7 α -type and 7 β -type subunits. Upon stimulation with IFN- γ , the three active β subunits, β 1 (Y), β 2 (Z) and β 5 (X) are exchanged to their immunocounterparts, namely β 1i (LMP-2), β 2i (MECL-1) and β 5i (LMP-7). This results in a change in proteasomal substrate specificity. Some groups have reported increases in chymotrypsin-like and trypsin-like activities while others have reported no changes or even decreased activities after such treatment.

Citations: 7

[View Online »](#)

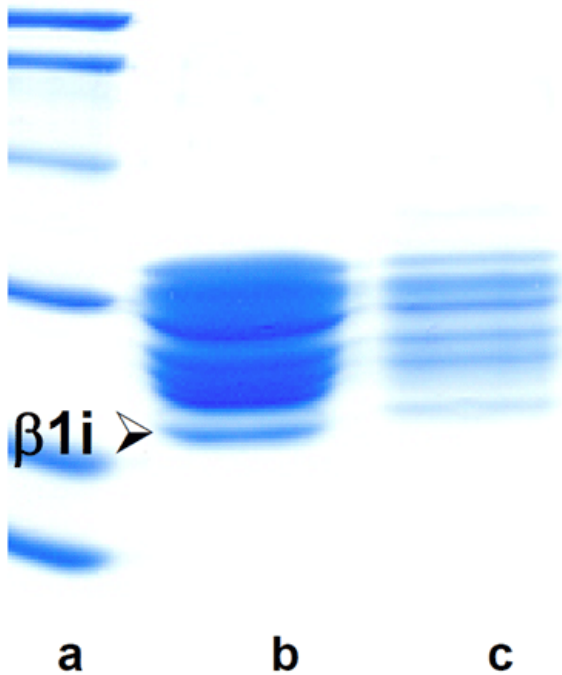
Ordering Information

[Order Online »](#)

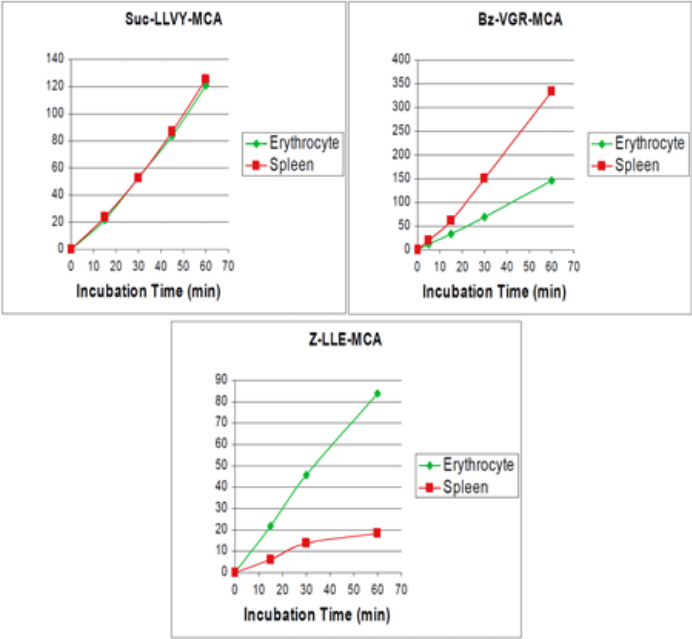
BML-PW9645-0050	50 μ g
-----------------	------------

Manuals, SDS & CofA

[View Online »](#)

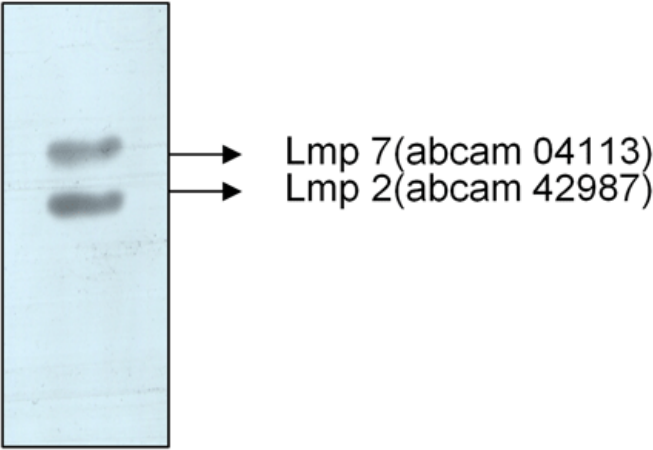


Coomassie stained gel after SDS-PAGE (12.5% acrylamide). Lane (a): molecular weight markers; lane (b) 20S immunoproteasome, spleen-derived (prod.no. BML-PW9645); lane (c): 20S proteasome, erythrocyte-derived (PW8720).



All activities were measured by use of 200μM substrate concentration (final concentration) all dissolved in TEAD buffer. Test system: 20μl enzyme solution (640ng); 20μl substrate solution (incubation at 37°C); 200μl stop solution (100mM sodium chloroacetate dissolved in 30mM Na-acetate, 70mM acetic acid, pH 4.3). Fluorescence was measured in a microplate fluorimeter at 355nm excitation and 460nm emission.

Immunoblot



Western blot analysis of Immunoproteasome 20S (human), (purified) (Prod. No. BML-PW9645) probed with anti-Syntaxin 3 antibody (ab4113) and anti-Proteasome 20S LMP2 antibody (ab42987).

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Once thawed the material can be stored at +4°C for up to 3 months, or long term at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Liquid. In TEAD buffer (20mM TRIS/HCl, 1mM sodium azide, 1mM DTT, pH 7.4). Contains 50% glycerol.
Purity Detail	Highly purified. All starting material has been tested and found to be negative for hepatitis B surface antigen, human immunodeficiency virus type 1 antigens, and antibodies against human immunodeficiency viruses type 1 and 2, and hepatitis C virus.
Source	Isolated from human spleen.

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)