

Granzyme B (human), (recombinant)

Granzyme B is a component of the cytolytic granules of cytotoxic T lymphocytes and natural killer cells.

Citations: 13

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE238-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Biological Activity	A serine protease with caspase-like substrate specificity.
Concentration	100 U/μl
Formulation	Liquid. In 50mM TRIS, pH 7.4, 100mM NaCl, 10% glycerol.
MW	25.4 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Source	Produced in yeast. Recombinant human granzyme B (aa 21-247).
Specific Activity	One unit hydrolyzes 1 pmol/min of Ac-IEPD-pNA (200 μM, Prod. No. BML-P133) at 30°C.
UniProt ID	P10144

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)