

Granzyme A (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-118-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -80°C. Do not dilute to more than 0.2mg/ml. Do not store dilute solutions (1-1000ng/ml) for more than 24 hours.
Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Concentration	0.2mg/ml
Formulation	Liquid. In 0.05M MES, pH 6.0, containing 10% sucrose and 0.5M sodium chloride. Contains no preservatives.
MW	~52kDa (homodimer)
Purity	≥98% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Active human granzyme A.
Specific Activity	Refer to the lot specific CofA for specific activity.
UniProt ID	P12544

Last modified: December 17, 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)