

Granzyme A (human)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-200-605-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months when stored at -80°C. 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

Formulation	Liquid. In ~0.15M sodium chloride.
MW	~52kDa (fully glycosylated homodimer)
Purity Detail	Purified by differential solubilization and cation exchange chromatography. Dialyzed against 0.15M sodium chloride and concentrated.
Source	Human cytotoxic granules of NK-92 cells.
Specific Activity	~9.2U/μg protein. One unit is defined as the amount of enzyme that hydrolyzes 1nmole of Z-Lys-SBzl per min. at 25°C in 0.05M TRIS, pH 8.0, containing 0.15M NaCl, 0.01% Triton X-100, 0.3mM DTNB.
UniProt ID	P12544



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)