

# Granzyme K (human), (recombinant)

Granzyme K is a trypsin-like serine protease located in the granules of natural killer (NK) cells and cytotoxic T lymphocytes (CTLs). It plays a crucial role in inducing apoptosis and also contributes to extracellular processes, such as activating endothelial cells and regulating immune responses. Elevated levels of Granzyme K have been associated with a range of diseases, including viral and bacterial infections.

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-117A-C010	10µg
ALX-201-117A-C5010	5x10µg

## Manuals, SDS & CofA

[View Online »](#)

- **High Specific Activity:** Granzyme K is provided with a defined specific activity, ensuring consistent and reliable enzymatic performance, with detailed activity data available in the lot-specific Certificate of Analysis (CoA).
- **Native Source and Activity:** Recombinant human granzyme K produced in *E. coli*, preserving native enzymatic function for accurate biological studies.
- **Broad Assay Compatibility:** Optimized to hydrolyze Z-Lys-SBzl in a range of standard assay conditions, facilitating diverse experimental applications.
- **Exceptional Purity:** Purity level of ≥98% (SDS-PAGE), ensuring minimal interference from contaminants in sensitive applications.

## Handling & Storage

Use/Stability	Dilute solutions (e. g. 1-1000ng/ml) should be used within 24 hours.
Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. 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 0.05 M sodium phosphate buffer, containing 0.5 M NaCl and 10% glycerol. Contains no preservatives.
MW	~26kDa
Purity	≥95% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Active human granzyme K.
Specific Activity	One unit is defined as the amount of enzyme that hydrolyzes 1nmol Z-Lys-SBzl per min. at 25°C in 0.05M TRIS, pH 8.0, containing 0.15M NaCl, 0.01% Triton X-100 and 0.3mM DTNB. Specific activity provided in lot specific CoA.
UniProt ID	P49863

Last modified: January 9, 2025



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)