

# Cathepsin G (human neutrophils), (purified)

Cathepsin G, a serine protease found in polymorphonuclear neutrophils (PMNs), functions in inflammation.

Citations: 11

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-SE283-0100	100µg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application Notes** Useful for inhibitor screening, kinetic studies, and cleavage of substrates *in vitro* or in tissue culture.

**CAS** 56645-49-9

**Formulation** Lyophilized.

**Purity** ≥95% by SDS-PAGE

**Reconstitution** Reconstitute with 50 mM sodium acetate, pH5.5, containing 150 mM NaCl.

**Source** Isolated from human neutrophils. Active enzyme. Negative for HbsAg, anti-HBc, anti-HCV, and anti-HIV-1 and -2.

**Specific Activity** 2 to 4 units per mg protein. One unit of activity is defined as the amount of enzyme that hydrolyzes one micromole of Suc-Ala-Ala-Pro-Phe-pNA (1mM) per minute at 25°C in 160mM TRIS-HCl, pH 7.4, containing 1.6M sodium chloride.

**UniProt ID** P08311



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-  
fr@enzolifesciences.com](mailto:info-fr@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)