

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

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-
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)