

Neutrophil elastase (human), (purified)

Major inflammatory protease released by neutrophils

Functions in inflammation. Can cleave a variety of proteins, including fibrous elastin, collagen, clotting factors, TNF- α , and α 1-antichymotrypsin. Because of its potential roles in inflammatory diseases such as emphysema and cystic fibrosis, it is a target for drug discovery.

Neutrophil elastase, a serine protease found in polymorphonuclear neutrophils (PMNs), functions in inflammation.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE284-0100	100 μ g
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. This enzyme is stable when stored under the above conditions. Procedures such as dilution of the enzyme followed by refreezing could lead to loss of activity.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Leukocyte elastase, LE
Application Notes	Useful for inhibitor screening, kinetic studies, and cleavage of substrates <i>in vitro</i> or in tissue culture.
Purity	≥95% (SDS-PAGE). Although this is in highly purified form, and the source material tested negative for HbsAg, anti-HBc, anti-HCV, and anti-HIV-1 and -2, handle this product with all due precautions.
Reconstitution	Reconstitute with 50mM sodium acetate, pH 5.5, with 150mM sodium chloride to the original volume, followed by gentle mixing to ensure adequate homogenization.
Source	Isolated from human neutrophils. Active enzyme.
Specific Activity	20-22 U/mg protein. One unit will hydrolyze 1.0 μmole of MeOSuc-Ala-Ala-Pro-Val-pNA (Prod. No. BML-P213) per minute, at 25°C in 100mM Tris-HCl, pH 7.5 with 500 mM NaCl.
UniProt ID	P08246



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)