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

100µg

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 in vitro 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

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com