

Neutrophil elastase colorimetric drug discovery kit

It is a complete assay system designed to screen inhibitors of neutrophil elastase, providing enough material to perform at least 96 assays. Neutrophil elastase, also known as leukocyte elastase (LE), is a serine protease found in polymorphonuclear neutrophils (PMNs) that functions in inflammation. It can cleave a variety of proteins, including fibrous elastin, collagen, clotting factors, TNF- α , and α 1-antichymotrypsin. It has been shown to be taken up by tumor cells and accelerate tumor growth through its degradation of insulin receptor substrate-1 (IRS-1). Because of its potential roles in cancer and inflammatory diseases such as emphysema, cystic fibrosis, and atherosclerosis, it is a target for drug discovery. The kit uses the chromogenic substrate MeOSuc-Ala-Ala-Pro-Val-pNA; cleavage of p-nitroaniline (pNA) from the substrate increases absorbance at 405nm. A neutrophil elastase inhibitor, elastatinal, is included for use as a control.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK497-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

Figure 3. Percent inhibition of HNE by Elastatinal

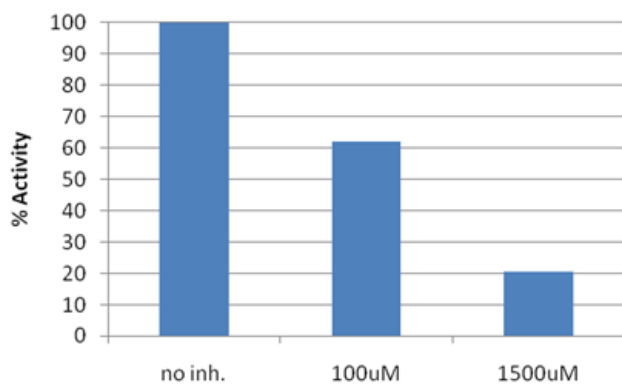


Figure 2. Inhibition of HNE by Elastatinal

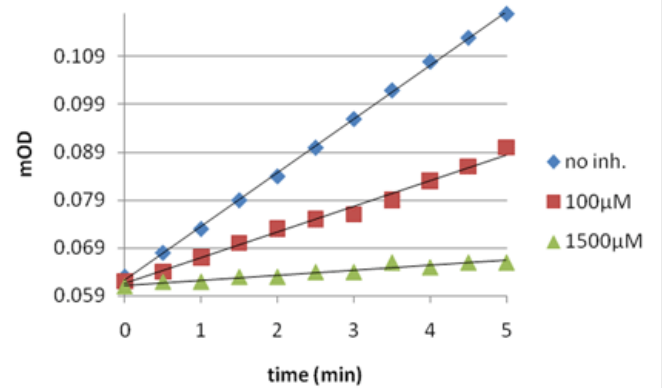
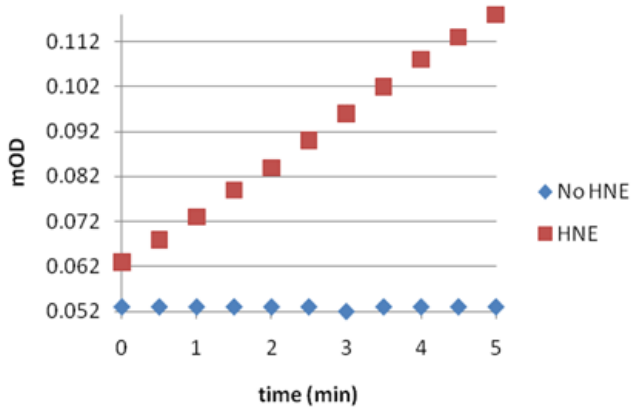


Figure 1. Plot of OD vs. time



Handling & Storage

Use/Stability	Substrates and inhibitor are in DMSO and are stable for up to 6 months. Neutrophil elastase activity decreases with time, even at -80°C.
Handling	Keep frozen at all times.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Activity assay, Colorimetric detection, HTS
Application Notes	Screening of neutrophil elastase inhibitors.
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	<p>Neutrophil elastase (human), (purified) (Prod. No. BML-SE284-9090)</p> <p>pNA substrate (MeOSuc-AAPV-pNA) (Prod. No. BML-P213-9090)(60µl; 20mM in DMSO)</p> <p>pNA calibration standard (p-nitroaniline) (Prod. No. BML-KI106)(1ml; 50µM in assay buffer)</p> <p>Elastatinal (inhibitor) (Prod. No. BML-PI103-9090)(10µl; 20mM in DMSO)</p> <p>Neutrophil elastase assay buffer (Prod. No. BML-KI581)(20ml; 100mM HEPES, pH 7.25, 500mM NaCl, 0.05% Tween-20; liquid in a screw-cap plastic bottle)</p> <p>96-well microplate (Prod. No. BML-KI101)</p>
UniProt ID	P08246

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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