

Neutrophil Elastase Activity Assay (fluorometric)

A plate based fluorometric assay to measure elastase activity in a variety of biological sample types with a sensitivity below 50 μ U in a sample well.

Neutrophil elastase is a key component of the immune response. The primary function of neutrophil elastase is to degrade bacterial components at the site of infection, contributing to the breakdown of the bacterial cell wall. In inflammatory disease conditions, neutrophil elastase becomes dysregulated, being released inappropriately by the neutrophils and degrading ECM proteins which are essential for tissue integrity and function. This is seen in conditions such as COPD, ARDS and autoimmune disorders like rheumatoid arthritis. The ability to selectively target neutrophil elastase with inhibitors in several disease states could mitigate excessive tissue destruction while preserving its activity in infections. The ability to measure neutrophil elastase activity is essential in this work. AkrisBio's Neutrophil Elastase Activity Assay is a simple, sensitive means of assessing elastase activity in a variety of biological sample types with a sensitivity below 50 μ U in a sample well.

Ordering Information

[Order Online »](#)

AKR-MA-0156	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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-be@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