

NETosis Assay Kit ab235979

★★★★★ [1 Abreviews](#) [2 Images](#)

Overview

Product name	NETosis Assay Kit
Detection method	Colorimetric
Sample type	Suspension cells
Assay type	Quantitative
Range	0.56 mU/well - 36 mU/well
Species reactivity	Reacts with: Human
Product overview	<p>NETosis Assay Kit (ab235979) provides a simple and fast method for studying the process of NETosis <i>ex vivo</i>.</p> <p>In this kit, primary neutrophils are stimulated to release NETs with either PMA or a calcium ionophore (both included). Unbound neutrophil elastase is washed away following NET generation. Following digest of NET DNA by S7 nuclease, the supernatant containing neutrophil elastase is added to a substrate, which is selectively cleaved by elastase to yield a 4-nitroaniline product that adsorbs light at 405 nm. Enough reagents are provided to test 24 sample conditions for NET production, with analysis in duplicate.</p> <p>Note: This kit does not depend upon the DNA component of neutrophil extracellular traps (NETs), as DNA release can occur independently of NETosis.</p>
Platform	Microplate reader

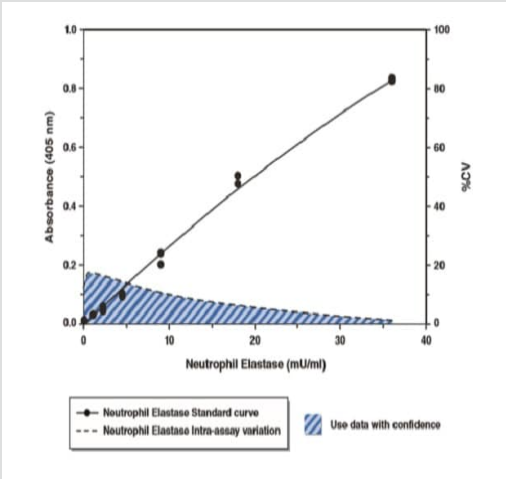
Properties

Storage instructions Please refer to protocols.

Components	1 kit	1 kit
96-Well Plate (Colorimetric Assay)	1 unit	1 unit
96-Well Plate Cover	1 unit	1 unit
A-23187 (25 mM) Assay Reagent	1 x 50µl	1 x 50µl

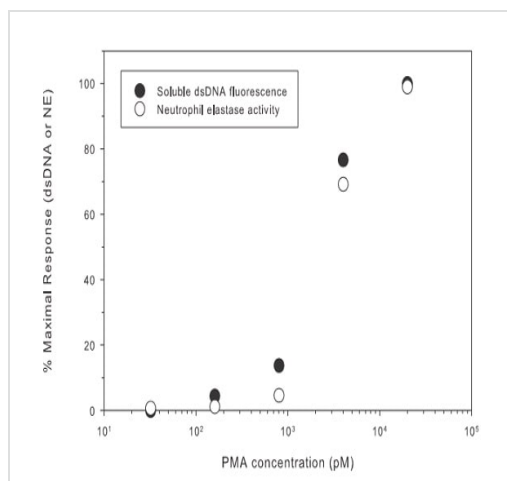
Components	1 kit	1 kit
Bovine Serum Albumin Assay Reagent	1 x 5g	1 x 5g
Calcium Chloride (1 M) Assay Reagent	1 x 1ml	1 x 1ml
Cell-Based PMA (1 mM)	1 x 50µl	1 x 50µl
EDTA (500 mM) Assay Reagent	1 x 1ml	1 x 1ml
Human Neutrophil Elastase Assay Reagent	1 x 50µl	1 x 50µl
NET Assay Neutrophil Elastase Substrate	2 vials	2 vials
S7 Nuclease Assay Reagent	1 x 50µl	1 x 50µl

Images



Human neutrophil elastase standard curve.

Example standard curve



Measurement of released neutrophil elastase parallels measurement of released dsDNA.

Human neutrophils were treated with PMA for 4 hours, washed and treated with S7 Nuclease for 15 minutes. The supernatant from each well was sampled and assayed for neutrophil elastase according to the protocol booklet. The supernatant was also tested for the presence of soluble dsDNA using a green fluorescence probe.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors