

Human Neutrophil Elastase ELISA Kit ab270204

Recombinant SimpleStep ELISA<sup>®</sup>

1 References 6 Images

Overview

Product name Human Neutrophil Elastase ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Serum	4			4.1%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Serum	8			3.2%

Sample type Serum, Cell Lysate, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 15.8 pg/ml

Range 47 pg/ml - 3000 pg/ml

Recovery	Sample specific recovery		
	Sample type	Average %	Range
	Serum	112	105% - 120%
	Cell culture extracts	116	115% - 117%
	Cell culture media	106	103% - 109%
	Hep Plasma	108	105% - 113%
	EDTA Plasma	111	103% - 118%
	Cit plasma	102	100% - 105%

<b>Assay time</b>	1h 30m
<b>Assay duration</b>	One step assay
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	<p>Human Neutrophil Elastase ELISA Kit (ab270204) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Neutrophil Elastase protein in cell culture media, cell lysate, edta plasma, serum, cit plasma, and hep plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Neutrophil Elastase with 15.8 pg/ml sensitivity.</p> <p>SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:</p> <ul style="list-style-type: none"> <li>- Single-wash protocol reduces assay time to 90 minutes or less</li> <li>- High sensitivity, specificity and reproducibility from superior antibodies</li> <li>- Fully validated in biological samples</li> <li>- 96-wells plate breakable into 12 x 8 wells strips</li> </ul> <p>A 384-well SimpleStep ELISA® microplate (<a href="#">ab203359</a>) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.</p>
<b>Notes</b>	<p>Neutrophil Elastase, also known as bone marrow serine protease, Elastase-2, human leukocyte Elastase, and medullasin is a serine protease in the peptidase S1 family and is encoded by the <i>ELANE</i> gene. Neutrophil Elastase is expressed as a zymogen by neutrophils and macrophages during inflammation, where it breaks down bacterial and host membrane proteins. Neutrophil Elastase can localize to Neutrophil extracellular traps (NETs) by the protein's affinity for DNA.</p> <p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.</p> <p>It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
<b>Platform</b>	Pre-coated microplate (12 x 8 well strips)

## Properties

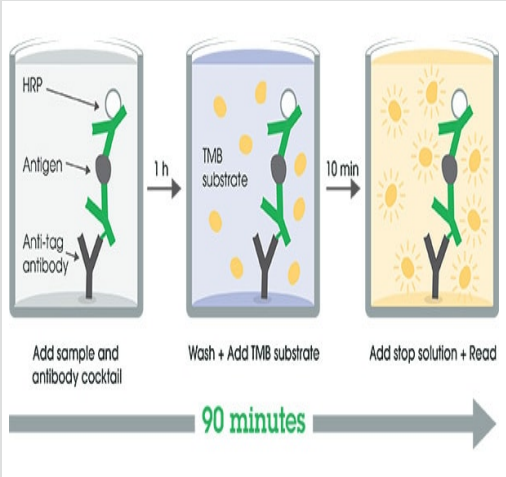
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Human Neutrophil Elastase Capture Antibody	1 x 600µl
10X Human Neutrophil Elastase Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml

Components	1 x 96 tests
Antibody Diluent 4BI	1 x 6ml
Human Neutrophil Elastase Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent 25BS	1 x 20ml
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

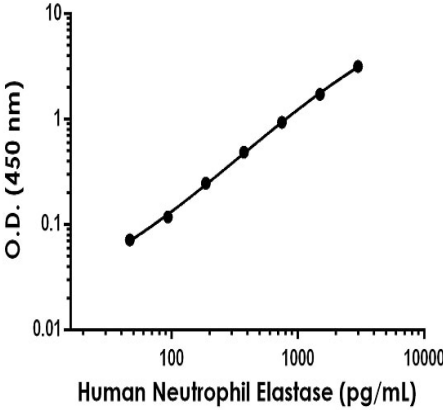
<b>Function</b>	Modifies the functions of natural killer cells, monocytes and granulocytes. Inhibits C5a-dependent neutrophil enzyme release and chemotaxis.
<b>Tissue specificity</b>	Bone marrow cells.
<b>Involvement in disease</b>	<p>Defects in ELANE are a cause of cyclic haematopoiesis (CH) [MIM:162800]; also known as cyclic neutropenia. CH is an autosomal dominant disease in which blood-cell production from the bone marrow oscillates with 21-day periodicity. Circulating neutrophils vary between almost normal numbers and zero. During intervals of neutropenia, affected individuals are at risk for opportunistic infection. Monocytes, platelets, lymphocytes and reticulocytes also cycle with the same frequency.</p> <p>Defects in ELANE are the cause of neutropenia severe congenital autosomal dominant type 1 (SCN1) [MIM:202700]. SCN1 is a disorder of hematopoiesis characterized by a maturation arrest of granulopoiesis at the level of promyelocytes with peripheral blood absolute neutrophil counts below <math>0.5 \times 10^9/l</math> and early onset of severe bacterial infections.</p>
<b>Sequence similarities</b>	<p>Belongs to the peptidase S1 family. Elastase subfamily.</p> <p>Contains 1 peptidase S1 domain.</p>

## Images



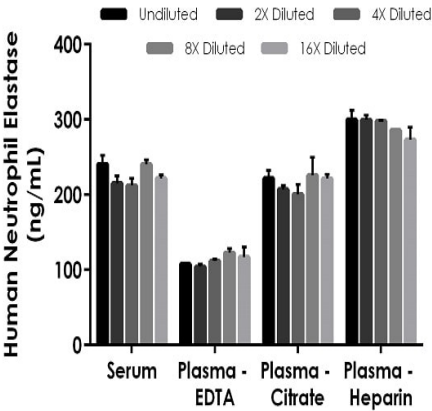
SimpleStep ELISA Protocol Diagram

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



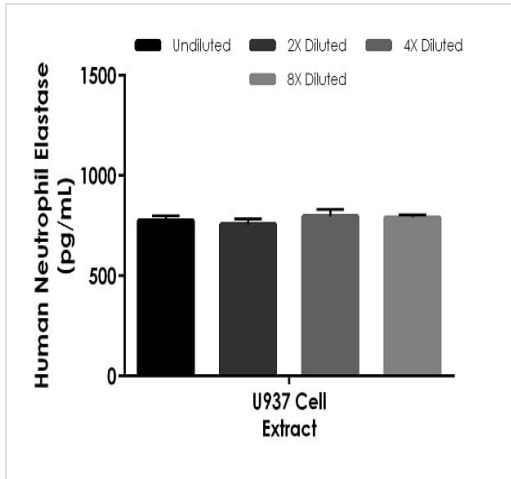
The Neutrophil Elastase standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean  $\pm$  SD) are graphed.

Example of human Neutrophil Elastase standard curve in Sample Diluent 1X Cell Extraction Buffer PTR.



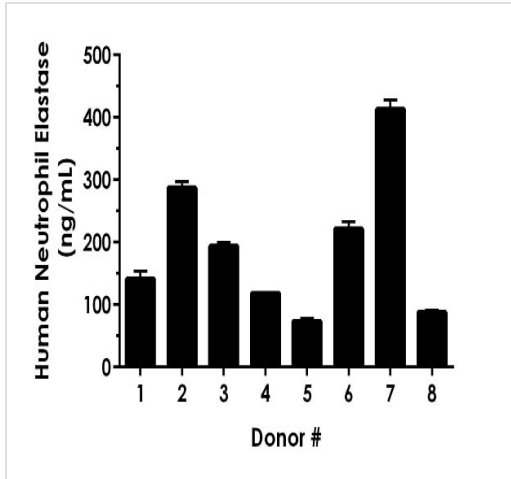
The concentrations of Neutrophil Elastase were measured in duplicates, interpolated from the target standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1:100, plasma (EDTA) 1:50, Plasma (citrate) 1:100, and plasma (heparin) 1:200. The interpolated dilution factor corrected values are plotted (mean  $\pm$  SD,  $n=2$ ). The mean target concentration was determined to be 225 ng/mL in serum, 113 ng/mL in plasma (EDTA), 215 ng/mL in plasma (citrate), and 291 ng/mL in plasma (heparin).

Interpolated concentrations of native Neutrophil Elastase in human serum, plasma (EDTA), plasma (citrate), plasma (heparin).



Interpolated concentrations of native Neutrophil Elastase in human U937 cell extract based on a 2.5 µg/mL extract load.

The concentrations of Neutrophil Elastase were measured in duplicate and interpolated from the Neutrophil Elastase standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean  $\pm$  SD, n=2).



Serum from eight individual healthy Neutrophil Elastase gender donors was measured in duplicate.

Interpolated dilution factor corrected values are plotted (mean  $\pm$  SD, n=2). The mean Neutrophil Elastase concentration was determined to be 192 ng/mL with a range of 74 – 413 ng/mL.

Powered by  
recombinant antibodies



**Research with confidence**  
Consistent and reproducible results



**Long-term and scalable supply**  
Recombinant technology



**Success from the first experiment**  
Confirmed specificity



**Ethical standards compliant**  
Animal-free production

Sandwich ELISA - Human Neutrophil Elastase  
ELISA Kit (ab270204)

To learn more about the advantages of recombinant antibodies see [here](#).

**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