

ab325362 – Human Neutrophil Elastase SimpleStep ELISA® Kit, Chemiluminescent

For the quantitative measurement of Neutrophil Elastase in human serum, plasma (heparin), plasma (EDTA), plasma (citrate), cell culture supernatant, CSF, and cell extract.

For research use only - not intended for diagnostic use.

Patent Pending

For overview, typical data and additional information please visit: www.abcam.com/ab325362

Storage and Stability: Store kit at 2-8°C immediately upon receipt. Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Standard Preparation and Reagent preparation sections.

Limitations: Typical standard curve, sensitivity, CSF linearity and recovery, and disease-state sample data were determined using Chemiluminescent detection. Unless stated, all remaining data was collected using the colorimetric version of this kit (ab270204).

Materials Supplied

Item	Quantity	Storage Condition
Human Neutrophil Elastase Capture Antibody 10X	600 µL	+4°C
Human Neutrophil Elastase Detector Antibody 10X	600 µL	+4°C
Human Neutrophil Elastase Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 4BI	6 mL	+4°C
Cell Extraction Buffer PTR 5X	10 mL	+4°C
Sample Diluent 25BS	20 mL	+4°C
Sample Diluent NS	12 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
ChemiHRP Reagent A	3 mL	+4°C
ChemiHRP Reagent B	3 mL	+4°C
SimpleStep Pre-Coated Black 96-Well Microplate	96 Wells	+4°C
Plate Seal	1	+4°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Luminometer with the following settings: 0.5-1 second/well read time; summation mode (all wavelengths).
- Method for determining protein concentration (BCA assay recommended).
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Plate shaker for all incubation steps.
- Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.

Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

Sample Diluent 25BS may contain precipitate, this is normal. If precipitate is not dissolved by gentle mixing, the precipitate may be dissolved by gentle warming and mixing at 37°C for 10 minutes. If precipitate remains, gently spin down and avoid visible precipitates when pipetting.

1X Cell Extraction Buffer PTR (For cell and tissue extracts only): Prepare 1X Cell Extraction Buffer PTR by diluting Cell Extraction Buffer PTR 5X and 50X Cell Extraction Enhancer Solution to 1X with deionized water. To make 10 mL 1X Cell Extraction Buffer PTR combine 8 mL deionized water and 2 mL Cell Extraction Buffer PTR 5X. Mix thoroughly and gently. If required protease inhibitors can be added.

1X Wash Buffer PT: Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 4BI. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent 4BI. Mix thoroughly and gently.

Lumi HRP Development Solution: Just prior to use, prepare Lumi HRP Development Solution by mixing equal volume of the ChemiHRP Reagent A and the ChemiHRP Reagent B. To make 3 mL of the Lumi HRP Development Solution combine 1.5 mL of ChemiHRP Reagent A and 1.5 mL of ChemiHRP Reagent B. Mix thoroughly and gently by inversion or slow pipetting (Avoid shaking or vortexing). Protect the prepared solution from light until use.

Standard Preparation

Always prepare a fresh set of standards for every use. Discard working standard dilutions after use as they do not store well. The following section describes the preparation of a standard curve for duplicate measurements (recommended).

- IMPORTANT:** If the protein standard vial has a volume identified on the label, reconstitute the Neutrophil Elastase protein standard by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the Neutrophil Elastase standard by adding 500 µL Diluent. Hold at room temperature for 10 minutes and mix gently. This is the 32,000 pg/mL Stock Standard Solution.
For serum and plasma samples measurements, reconstitute the Neutrophil Elastase protein standard by adding 25BS. For **cell culture supernatant and CSF** samples measurements, reconstitute the Neutrophil Elastase protein standard by adding Sample Diluent NS. For **cell extracts** samples measurements, reconstitute the Neutrophil Elastase protein standard by adding 1X Cell Extraction Buffer PTR.
- Label nine tubes, Standards 1–9.
- Use the same Sample Diluent as used to resuspend the Stock Standard to prepare the standard curve. Add 220 µL of Sample Diluent into tube number 1 and 150 µL of Sample Diluent into numbers 2-9.
- Use the **Stock Standard** to prepare the following dilution series. Standard #9 contains no protein and is the Blank control:

Standard #	Dilution Sample	Volume to Dilute (µL)	Volume of Diluent (µL)	Starting Conc. (pg/mL)	Final Conc. (pg/mL)
1	Stock Standard	100	220	32,000	10,000
2	Standard#1	75	150	10,000	3,333.3
3	Standard#2	75	150	3,333.3	1,111.1
4	Standard#3	75	150	1,111.1	370.4
5	Standard#4	75	150	370.4	123.5
6	Standard#5	75	150	123.5	41.2
7	Standard#6	75	150	41.2	13.7
8	Standard#7	75	150	13.7	4.6
9	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	1:3,225 - 1:100
Plasma – Citrate	1:3,225 - 1:100
Plasma – EDTA	1:1,667 - 1:50
Plasma – Heparin	1:3,225 - 1:200
Cell Culture Media*	≤ 50%
Cerebrospinal Fluid	6.25 - 50%
U937 Cell Extract	0.31 - 2.5 µg/mL

*Based on spiked sample

Serum Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Dilute samples at least 1:100 into Sample Diluent 25BS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

Plasma Collect plasma using citrate, EDTA, or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Dilute plasma Citrate samples at least 1:100, plasma EDTA samples at least 1:50, and plasma citrate samples at least 1:200 into Sample Diluent 25BS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

Cell Culture Supernatants Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Dilute samples at least 1:2 into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Cerebrospinal Fluid (CSF) Dilute cerebrospinal fluid at least 1:2 into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Preparation of extracts from cell pellets Collect non-adherent cells by centrifugation or scrape to collect adherent cells from the culture flask. Typical centrifugation conditions for cells are 500 x g for 5 minutes at 4°C. Rinse cells twice with PBS. Solubilize pellet at 2×10^7 cell/mL in chilled 1X Cell Extraction Buffer PTR. Incubate on ice for 20 minutes. Centrifuge at 18,000 x g for 20 minutes at 4°C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay. Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

Preparation of extracts from adherent cells by direct lysis (alternative protocol) Remove growth media and rinse adherent cells 2 times in PBS. Solubilize the cells by addition of chilled 1X Cell Extraction Buffer PTR directly to the plate (use 750 µL - 1.5 mL 1X Cell Extraction Buffer PTR per confluent 15 cm diameter plate). Scrape the cells into a microfuge tube and incubate the lysate on ice for 15 minutes. Centrifuge at 18,000 x g for 20 minutes at 4°C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay. Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

Plate Preparation

The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.

Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.

For each assay performed, a minimum of two wells must be used as the zero control.

For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).

Differences in well absorbance or "edge effects" have not been observed with this assay.

Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use.

We recommend that you assay all standards, controls and samples in duplicate

1. Prepare all reagents, working standards, and samples as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
3. Add 50 µL of all sample or standard to appropriate wells.
4. Add 50 µL of the Antibody Cocktail to each well.
5. Seal the plate and incubate for 30 minutes at room temperature on a plate shaker set to 750 rpm.
6. Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 30 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
7. Add 50 µL of prepared Lumi HRP Development Solution to each well and incubate for 1 minute in the dark on a plate shaker set to 750 rpm. Further optimization of incubation time vs signal strength can be performed if needed. Avoid introducing bubbles into the wells.
8. Measure the produced light of each well using a microplate luminometer with the following settings: 0.5-1 second/well read time in summation mode (all wavelengths). Relative light unit (RLU) readings may vary between luminometer models. It is recommended to configure instrument settings according to the manufacturer's specifications. Note: Relative light unit (RLU) values may change over the course of the 15-minute reading window.
9. Analyze the data as described below.

Mode:	Luminescence
Instrument settings:	Endpoint
Detection Mode:	All wavelengths
Read Time:	0.5-1 sec
Read:	Top

Note For microplate readers with Pre-Read Optimization option, the Read Height as well as Microplate Optimization is recommended before the first read.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

Copyright © 2025 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.

Version 1b | 2025-12-11

ab325362 – Human Neutrophil Elastase SimpleStep ELISA® Kit, Chemiluminescent

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of human Neutrophil Elastase.

The standard protein in this kit is mature full-length human Neutrophil Elastase.

Native signal was detected in serum, plasma (heparin), plasma (EDTA), plasma (citrate), CSF, and cell extract.

Spiked protein experiments were used to validate cell culture supernatant sample types.

Saliva, urine, milk, and tissue extract samples have not been tested with this kit.

SPECIES REACTIVITY

Other species reactivity was determined by measuring 1:500 serum samples of various species, interpolating the protein concentrations from the human standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in human serum assayed at the same dilution.

No signal was observed for the following species: Mouse, Rat, Cow

Other species reactivity not determined.

CALCULATION

- Preconfigured protocols are available when using SoftMax Pro software from Molecular Devices.
- Calculate the average chemiluminescence value for the blank control (zero) standards. Subtract the average blank control standard chemiluminescence value from all other chemiluminescence values.
- Create a standard curve by plotting the average blank control subtracted chemiluminescence value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.
- Note: Most chemiluminescence reader software or graphing software will plot these values and fit a curve to the data. A four-parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4-parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.
- Determine the concentration of the target protein in the sample by interpolating the blank control subtracted chemiluminescence values against the standard curve. Multiply the resulting value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.
- Samples generating chemiluminescence values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at chemiluminescence values less than that of the lowest standard should be retested in a less dilute form.

TYPICAL DATA

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed

Standard Curve Measurements			
Concentration (pg/mL)	RLU		Mean RLU
	1	2	
0	2,601	2,601	2,601
4.6	3,588	3,389	3,489
13.7	5,810	5,856	5,833
41.2	13,011	13,463	13,237
123.5	36,188	35,979	36,084
370.4	111,490	113,220	112,355
1,111.1	334,630	340,850	337,740
3,333.3	969,140	946,980	958,060
10,000	3,098,400	2,961,900	3,030,150

Table 1. Example of human Neutrophil Elastase standard curve in Sample Diluent 25BS. The Neutrophil Elastase standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

Standard Curve Measurements			
Concentration (pg/mL)	RLU		Mean RLU
	1	2	
0	1,917	1,917	1,917
4.6	3,315	3,108	3,212
13.7	5,501	5,294	5,398
41.2	13,017	13,181	13,099
123.5	37,262	35,679	36,471
370.4	115,200	112,150	113,675
1,111.1	372,020	362,850	367,435
3,333.3	1,085,800	1,082,100	1,083,950
10,000	3,111,100	3,197,400	3,154,250

Table 2. Example of human Neutrophil Elastase standard curve in Sample Diluent NS. The Neutrophil Elastase standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

Standard Curve Measurements			
Concentration (pg/mL)	RLU		Mean RLU
	1	2	
0	3,202	3,202	3,202
4.6	4,604	4,449	4,527
13.7	6,750	7,121	6,936
41.2	13,958	13,199	13,579
123.5	37,208	34,896	36,052
370.4	110,640	107,800	109,220
1,111.1	327,770	337,820	332,795
3,333.3	1,058,100	1,058,300	1,058,200
10,000	3,164,200	3,186,500	3,175,350

Table 3. Example of human Neutrophil Elastase standard curve in 1X Cell Extraction Buffer PTR. The Neutrophil Elastase standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

TYPICAL SAMPLE VALUES

Sensitivity:

The minimal detectable dose (MDD) was determined by calculating the mean of zero standard replicates and adding 2 standard deviations then extrapolating the corresponding concentration.

Sample Diluent Buffer	N=	Minimal Detectable Dose
Sample Diluent 25BS	8	4.5 pg/mL
Sample Diluent NS	8	2.0 pg/mL
1X Cell Extraction Buffer PTR	16	4.2 pg/mL

Recovery

Three concentrations of Neutrophil Elastase were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
1:400 Serum	112	105 - 120
1:667 Plasma (citrate)	102	100 - 105
1:667 Plasma (EDTA)	111	103 - 118
1:667 Plasma (heparin)	108	105 - 113
50% Cell Culture Media*	106	103 - 109
12.5% Cerebrospinal Fluid	97	91 - 106
1 µg/mL U937 Cell Extract	116	115 - 117

*Media is RPMI 1640 containing 10% fetal calf serum.

Linearity of Dilution

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Native Neutrophil Elastase was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent 25BS.

Dilution Factor	Interpolated value	1:100 Human Serum	1:100 Human Plasma (Citrate)	1:50 Human Plasma (EDTA)	1:200 Human Plasma (Heparin)
Undiluted	pg/mL	2,409	2,217	2,159	1,499
	% Expected value	100	100	100	100
2	pg/mL	1,075	1,035	1,043	749
	% Expected value	89	93	97	100
4	pg/mL	530	501	560	373
	% Expected value	88	90	104	99
8	pg/mL	301	282	307	179
	% Expected value	100	102	114	95
16	pg/mL	139	139	147	85
	% Expected value	92	100	109	91

Neutrophil Elastase was measured in the following biological samples in a 2-fold dilution series. Native signal was measured in CSF and recombinant Neutrophil Elastase standard was spiked into cell culture media. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	50% Human CSF	50% Cell Culture Media*
Undiluted	pg/mL	49	4,988
	% Expected value	100	100
2	pg/mL	26	1,933
	% Expected value	106	79
4	pg/mL	14	1,326
	% Expected value	111	106
8	pg/mL	7	638
	% Expected value	107	102
16	pg/mL	ND	374
	% Expected value	ND	120

ND – Not Detected – below product dynamic range

*Based on spiked sample

Native Neutrophil Elastase was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in 1X Cell Extraction Buffer PTR.

Dilution Factor	Interpolated value	2.5 µg/mL U937 Cell Extract
Undiluted	pg/mL	775
	% Expected value	100
2	pg/mL	378
	% Expected value	98
4	pg/mL	199
	% Expected value	103
8	pg/mL	99
	% Expected value	102

Precision

Mean coefficient of variations of interpolated values of Neutrophil Elastase from a single concentration of serum within the working range of the assay.

	Intra-assay	Inter-assay
N=	4	8
CV (%)	4.1	3.2

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

Copyright © 2025 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.
Version 1b | 2025-12-11