

# Neutrophil Elastase Activity Assay Kit (Fluorometric) ab204730

★★★★☆ [1 Abreviews](#) [14 References](#) [2 图像](#)

### 概述

<b>产品名称</b>	Neutrophil Elastase Activity Assay试剂盒(Fluorometric)
<b>检测方法</b>	Fluorescent
<b>样品类型</b>	Plasma, Whole Blood, Purified protein
<b>检测类型</b>	Enzyme activity
<b>灵敏度</b>	1 ng
<b>检测时间</b>	0h 20m
<b>种属反应性</b>	<b>与反应:</b> Mammals, Other species
<b>产品概述</b>	<p>Neutrophil Elastase Activity Assay Kit (Fluorometric) ab204730 is based on the ability of neutrophil elastase (NE) to proteolytically cleave a synthetic substrate in order to release a fluorophore, AFC (Ex/Em 380/500 nm), which can be easily quantified by fluorescence microplate readers.</p> <p>This assay kit is simple, rapid and can detect as low as 1 ng of Neutrophil Elastase in a variety of samples.</p> <p>Neutrophil elastase assay protocol summary:</p> <ul style="list-style-type: none"> <li>- add samples and standards to wells</li> <li>- add reaction mix</li> <li>- analyze with a microplate reader every 2-3 min for 10-20 min</li> </ul>
<b>说明</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K383 Neutrophil Elastase Activity Assay kit (Fluorometric). K383-100 is the same size as the 100 test size of ab204730.</p> <p>Neutrophil Elastase (NE, EC 3.4.21.37, leukocyte elastase, ELANE, ELA2, elastase 2, neutrophil elaszym, or serine elastase) is a cytotoxic serine protease. It is stored in the azurophil granules of neutrophil granulocyte and is released following cell stimulation, e.g. by pathogens, immune complexes or chemotactic agents (PMA). When the extracellular NE concentration exceeds the buffering capacity of endogenous inhibitors, it causes degradation of a wide range of extracellular matrix proteins, including fibronectin, laminin, proteoglycans, collagens, and elastin.</p>
<b>平台</b>	Microplate reader

### 性能

## 存放说明

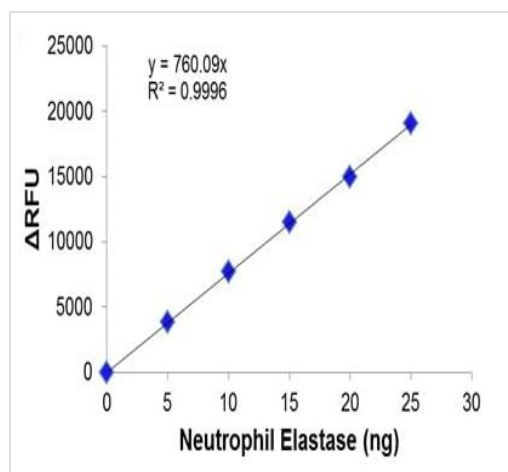
Store at -20°C. Please refer to protocols.

组件	100 tests
NE Assay Buffer	1 x 15ml
NE Dilution Buffer	1 x 1ml
Elastase Enzyme	1 x 1µg
Substrate VI	1 x 200µl

## 相关性

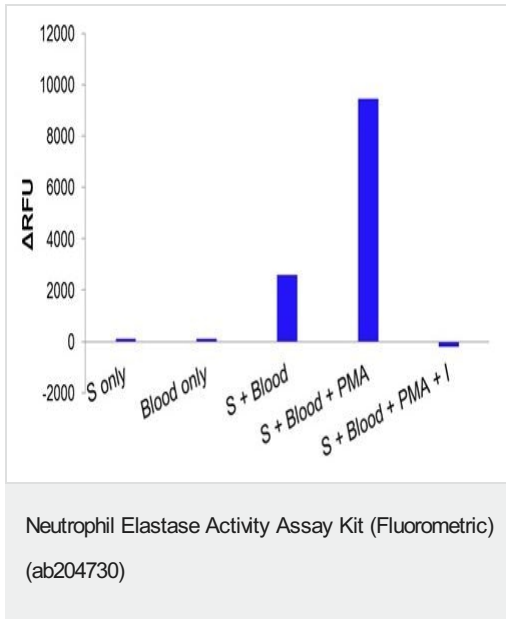
Neutrophil Elastase (NE) is an aggressive and cytotoxic 29 kDa serine protease stored mainly in the azurophil granules of neutrophil granulocytes. It plays a role in the degradation of a wide range of extracellular matrix proteins, including fibronectin, laminin, proteoglycans, collagens, and elastin. When the extracellular NE concentration exceeds the buffering capacity of endogenous inhibitors, it becomes implicated in the signs, symptoms and disease progression of inflammatory lung disorders via its role in the inflammatory process, mucus overproduction and lung tissue damage.

## 图片



Typical Neutrophil Elastase standard curve.

Neutrophil Elastase Activity Assay Kit (Fluorometric)  
(ab204730)



Neutrophil Elastase activity was measured in human blood, following removal of RBCs and stimulation of neutrophils by 100 nM PMA. S = Substrate, I = Inhibitor (60  $\mu$ M SPCK).

**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.cn/abpromise> or contact our technical team.

### Terms and conditions

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