

# Neutrophil Elastase Inhibitor Assay Kit ab118971

[13 References](#) [4 图像](#)

### 概述

<b>产品名称</b>	Neutrophil Elastase抑制剂Assay 试剂盒
<b>检测方法</b>	Fluorescent
<b>检测类型</b>	Quantitative
<b>检测时间</b>	0h 40m
<b>产品概述</b>	<p>In Abcam's Neutrophil Elastase Inhibitor Screening Kit, NE hydrolyzes a specific fluorescent substrate to release the fluorescent group, which can be detected at Ex/Em = 400/505 nm. In presence of a potent Neutrophil Elastase inhibitor, the hydrolyzation of substrate will be inhibited or stopped. The kit provides a rapid, simple, sensitive, and reliable test suitable as a high throughput screening assay of Neutrophil Elastase inhibitors. For comparison of the relative efficacy of test inhibitors, a control inhibitor, SPCK (Ki = 10 <math>\mu</math>M for human leukocyte elastase) is included.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
<b>说明</b>	This product is manufactured by BioVision, an Abcam company and was previously called K782 Neutrophil Elastase Inhibitor Screening Kit (Fluorometric). K782-100 is the same size as the 100 test size of ab118971.
<b>平台</b>	Microplate reader

### 性能

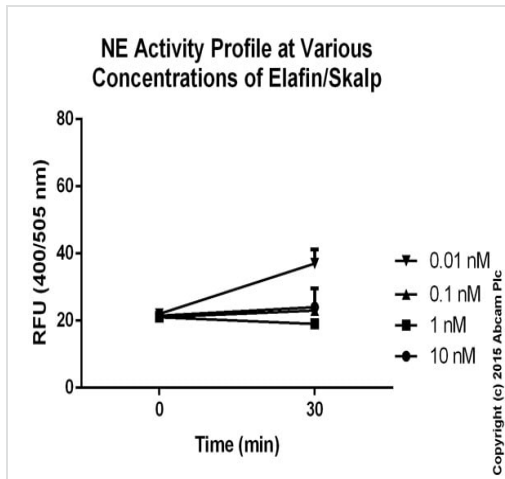
**存放说明** Store at -20°C. Please refer to protocols.

组件	100 tests
Assay Buffer XL	1 x 25ml
Elastase Inhibitor	1 x 100 $\mu$ l
Elastase Enzyme	1 vial
Substrate VI	1 x 0.2ml

**相关性** Neutrophil Elastase (NE) is an aggressive and cytotoxic 29 kDa serine protease stored mainly in

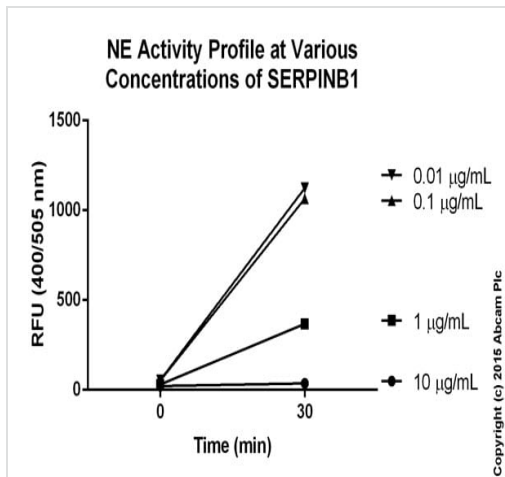
the azurophil granules of neutrophil granulocytes. It plays a role in the degradation of a wide range of extracellular matrix proteins, including fibronectin, laminin, proteoglycans, collagens, and elastin. When the extracellular NE concentration exceeds the buffering capacity of endogenous inhibitors, 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



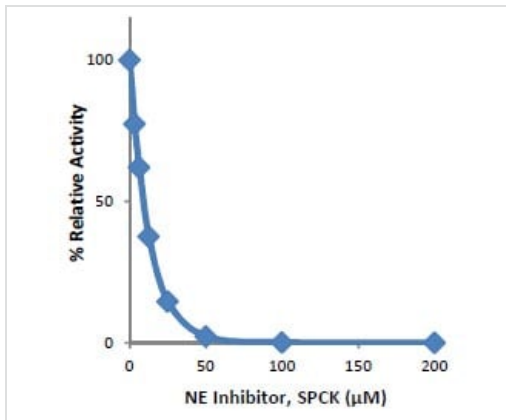
Neutrophil Elastase Activity Profile at Various Concentrations of Elafin/Skalp ([ab153768](#))

Neutrophil Elastase Inhibitor Screening Kit  
(ab118971)



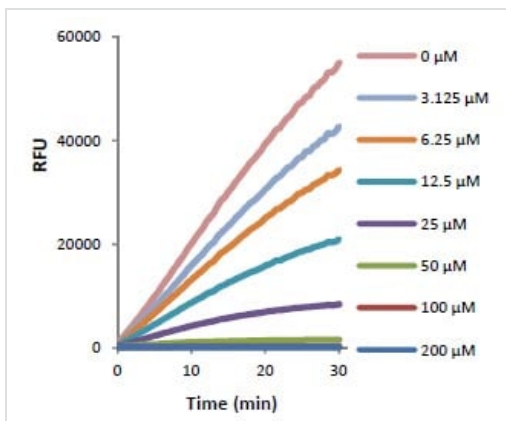
Neutrophil Elastase Activity Profile at Various Concentrations of SERPINB1 ([ab114774](#))

Neutrophil Elastase Inhibitor Screening Kit  
(ab118971)



Neutrophil Elastase Inhibition by SPCK

Functional Studies - Neutrophil Elastase Inhibitor Screening Kit (ab118971)



Neutrophil Elastase Activity Profile at Various Concentrations of SPCK

Functional Studies - Neutrophil Elastase Inhibitor Screening Kit (ab118971)

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

**Our Promise 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