

## Mouse eNOS ELISA Kit ab230938

**重组** SimpleStep ELISA

**2 References** [4 图像](#)

### 概述

**产品名称** 小鼠eNOS ELISA试剂盒

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
bEnd.3 cells	9			5%

批次间

样品	n	Mean	SD	CV%
bEnd.3 cells	3			5.04%

**样品类型** Cell culture extracts, Tissue Extracts, Mouse brain extract, Mouse muscle extract

**检测类型** Sandwich (quantitative)

**灵敏度** 5.58 pg/ml

**范围** 62.5 pg/ml - 4000 pg/ml

**回收率** 特定样本回收率

样品类型	平均%	范围
Cell culture extracts	102	101% - 103%
Tissue Extracts	96	95% - 98%
Mouse brain extract	96	95% - 97%
Mouse muscle extract	106	102% - 108%

**检测时间** 1h 30m

**实验步骤** One step assay

**种属反应性** 与反应: Mouse

## 产品概述

Mouse eNOS ELISA Kit (ab230938) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of eNOS protein in tissue extracts, mouse brain extract, mouse muscle extract, and cell culture extracts. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse eNOS with 5.58 pg/ml sensitivity.

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:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

## 说明

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. 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.

## 平台

Pre-coated microplate (12 x 8 well strips)

## 性能

### 存放说明

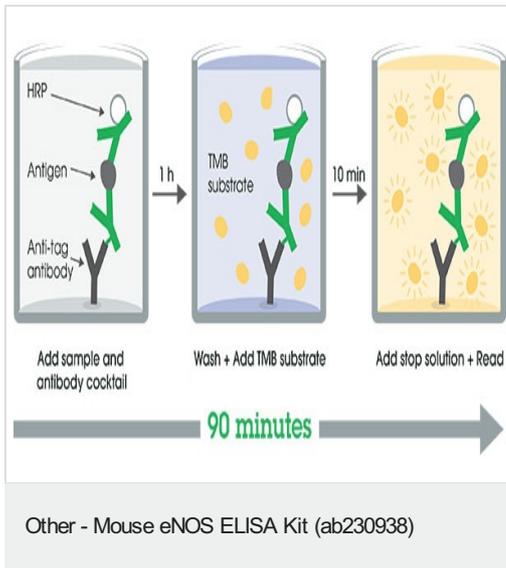
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Mouse eNOS Capture Antibody	1 x 600µl
10X Mouse eNOS Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent CPR	1 x 6ml
Mouse eNOS Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

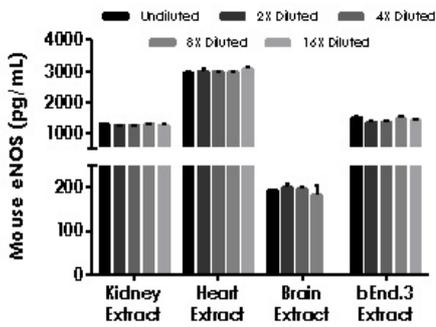
组件	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

<b>功能</b>	<p>Produces nitric oxide (NO) which is implicated in vascular smooth muscle relaxation through a cGMP-mediated signal transduction pathway. NO mediates vascular endothelial growth factor (VEGF)-induced angiogenesis in coronary vessels and promotes blood clotting through the activation of platelets.</p> <p>Isoform eNOS13C: Lacks eNOS activity, dominant-negative form that may down-regulate eNOS activity by forming heterodimers with isoform 1.</p>
<b>组织特异性</b>	Platelets, placenta, liver and kidney.
<b>疾病相关</b>	Variation in NOS3 seem to be associated with susceptibility to coronary spasm.
<b>序列相似性</b>	<p>Belongs to the NOS family.</p> <p>Contains 1 FAD-binding FR-type domain.</p> <p>Contains 1 flavodoxin-like domain.</p>
<b>翻译后修饰</b>	Phosphorylation by AMPK at Ser-1177 in the presence of Ca(2+)-calmodulin (CaM) activates activity. In absence of Ca(2+)-calmodulin, AMPK also phosphorylates Thr-495, resulting in inhibition of activity (By similarity). Phosphorylation of Ser-114 by CDK5 reduces activity.
<b>细胞定位</b>	Cell membrane. Membrane, caveola. Cytoplasm, cytoskeleton. Golgi apparatus. Specifically associates with actin cytoskeleton in the G2 phase of the cell cycle and which is favored by interaction with NOSIP and results in a reduced enzymatic activity.

## 图片

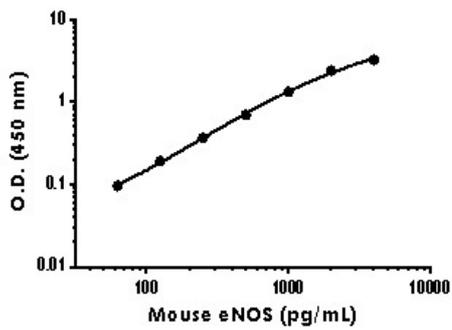


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Interpolated concentrations of native eNOS in mouse kidney extract, heart extract, brain extract, and bEnd.3 cell extract samples

The concentrations of eNOS were measured in duplicates, interpolated from the eNOS standard curves and corrected for sample dilution. Undiluted samples are as follows: kidney extract 1,000  $\mu\text{g/mL}$ , heart extract 1,000  $\mu\text{g/mL}$ , brain extract 250  $\mu\text{g/mL}$ , and bEnd.3 extract 25  $\mu\text{g/mL}$ . The interpolated dilution factor corrected values are plotted (mean  $\pm$  SD,  $n=2$ ). The mean eNOS concentration in the undiluted samples was determined to be 1,259  $\text{pg/mL}$  in kidney extract, 2,998  $\text{pg/mL}$  in heart extract, 194  $\text{pg/mL}$  in brain extract, and 1,419  $\text{pg/mL}$  in bEnd.3 cell extract.



Example of mouse eNOS standard curve in 1X Cell Extraction Buffer PTR

The eNOS 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.

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 - Mouse eNOS ELISA Kit  
(ab230938)

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.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