

### Nitric Oxide Assay Kit (Colorimetric) ab65328

★★★★★ [2 Abreviews](#) [99 References](#) [画像数 5](#)

医薬用外劇物

#### 製品の概要

製品名	Nitric Oxide Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate
アッセイタイプ	Quantitative
検出感度	1 nmol/well
検出範囲	1 nmol/well - 10 nmol/well
全工程の試験時間	1h 30m
種交差性	交差種: Mammals, Other species
製品の概要	Nitric Oxide Assay Kit (Colorimetric) ab65328 provides an accurate, convenient measure of total nitrate/nitrite in a simple two-step process.

In the nitric oxide assay protocol:

- the first step converts nitrate to nitrite by using nitrate reductase
  - the second step uses Griess Reagent to convert nitrite to a deep purple azo compound.
- The amount of the azochromophore accurately reflects nitric oxide amount in samples.

Nitric oxide assay protocol summary:

- add samples and standards to wells
- add nitrate reductase and enzyme cofactor and incubate for 1 hr at room temp to convert nitrate to nitrite
- add enhancer and incubate for 10 min at room temp
- add Griess Reagent R1 and Griess Reagent R2
- analyze with microplate reader

#### 特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K262 Nitric Oxide Colorimetric Assay Kit. K262-200 is the same size as the 2 96 tests size of ab65328.

If you are interested in a fluorometric detection kit, please check [Nitric Oxide Assay Kit \(Fluorometric\) \(ab65327\)](#).

For an assay kit to detect nitrite alone see [Griess Reagent Kit ab234044](#).

Nitric oxide (NO) plays an important role in neurotransmission, vascular regulation, immune response and apoptosis. NO is rapidly oxidized to nitrite and nitrate which are used to quantitate NO production.

## 製品の特性

## 保存方法

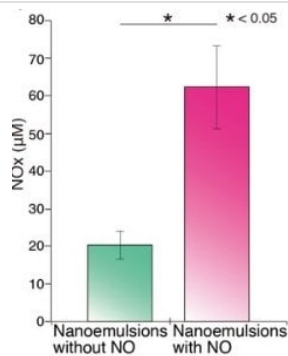
Store at -20°C. Please refer to protocols.

内容	2 x 96 tests
Assay Buffer XVII	1 x 30ml
Enhancer II	1 vial
Enzyme Cofactor	1 vial
Griess Reagent I	1 x 10ml
Griess Reagent II	1 x 10ml
96-Well Clear Plate	2 units
Nitrate Reductase I	1 vial
Nitrate Standard	1 vial
Nitrite Standard	1 vial
Microplate Sealing Film	2 units

## 関連性

Nitric oxide (NO) is a key vertebrate biological messenger, playing an important role in neurotransmission, vascular regulation, immune responses and apoptosis. NO, also known as "endothelium-derived relaxing factor" or "EDRF", is synthesized from L-arginine, oxygen and NADPH by various NO synthases. Most of the NO in the cell is oxidized to nitrite ( $\text{NO}_2^-$ ) and nitrate ( $\text{NO}_3^-$ ), and therefore the concentrations of these anions are generally as a quantitative measure of NO production.

## 画像



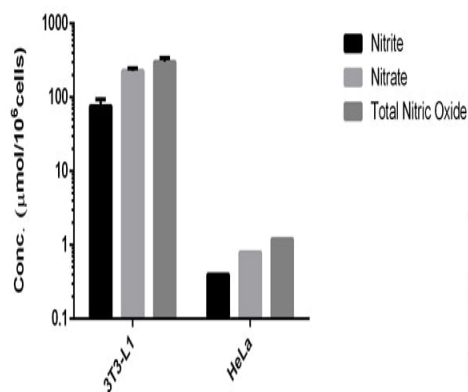
Nam, Eunryel et al.  
Scientific reports vol. 8,1  
7312. (2018)

#### Functional Studies - Nitric Oxide Assay Kit

(Colorimetric) (ab65328)

Nam, Eunryel et al., Scientific reports vol. 8,1 7312., Fig 4, doi:10.1038/s41598-018-25786-x

Blood NOx concentration measured with ab65328. Blood NOx concentration after application of the NE with NO. The level of blood NOx concentration increased 3-fold after the application of the NE with NO compared with control group.

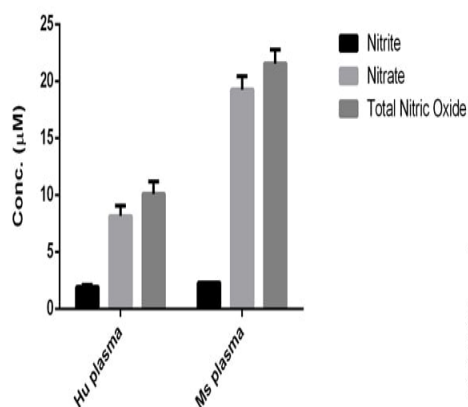


Copyright (c) 2014 Abcam plc

#### Functional Studies - Nitric Oxide Colorimetric Assay

Kit (ab65328)

Nitrite and nitrate measured in cell lysates showing quantity (micromol) per 10<sup>6</sup> cells

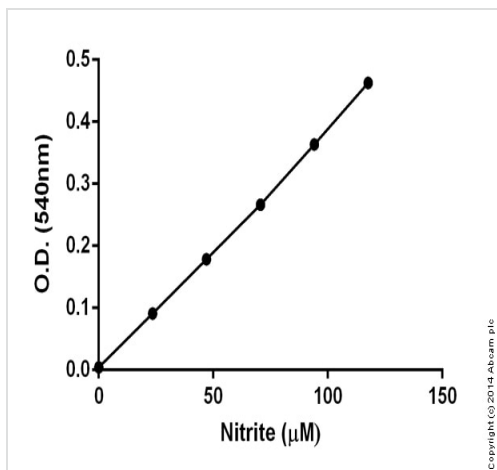


Copyright (c) 2014 Abcam plc

#### Functional Studies - Nitric Oxide Colorimetric Assay

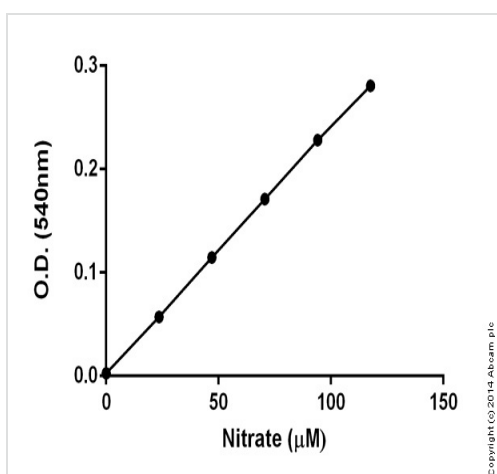
Kit (ab65328)

Nitrite and nitrate measured in biologicals showing concentration (micromolar)



Standard curve (colorimetric) : mean of duplicates (+/-SD) with background readings subtracted

Functional Studies - Nitric Oxide Colorimetric Assay  
Kit (ab65328)



Standard curve (Colorimetric) : mean of duplicates (+/-SD) with background readings subtracted

Functional Studies - Nitric Oxide Colorimetric Assay  
Kit (ab65328)

**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.co.jp/abpromise> or contact our technical team.

## **Terms and conditions**

---

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