

Product datasheet

Nitric Oxide Assay Kit (Fluorometric - Orange) ab219932

[2 References](#) [2 Images](#)

Overview

Product name	Nitric Oxide Assay Kit (Fluorometric - Orange)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Nitric Oxide Assay Kit (Fluorometric - Orange) (ab219932) is a sensitive fluorometric assay to monitor intracellular nitric oxide (NO) levels in live cells. The assay uses an orange dye that can react with NO to generate a bright orange fluorescent product that can be easily detected at Ex/Em = 540/590 nm, using the same filter set as Cy3[®] or TRITC.</p> <p>The assay can be detected by fluorescence microscopy, microplate fluorometry or high-content imaging. It can be easily adapted to use in 384-well microplate format.</p>
Notes	<p>Nitric oxide (NO) is an important biological regulator involved in numbers of physiological and pathological processes. Altered NO production is implicated in various immunological, cardiovascular, neurodegenerative and inflammatory diseases. As a free radical, NO is rapidly oxidized and there is relatively low concentrations of NO existing <i>in vivo</i>. It has been challenging to detect and understand the role of NO in biological systems.</p>
Platform	Microplate reader, Fluor. microscope, Flow cyt.

Properties

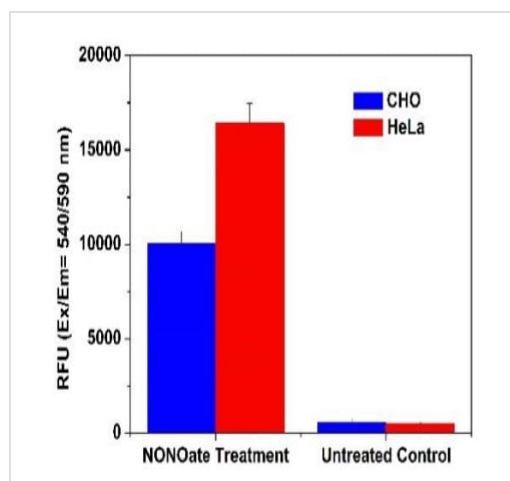
Storage instructions Store at -20°C. Please refer to protocols.

Components	200 tests
500X NO Orange Dye	1 x 50µl
Assay Buffer I	1 x 20ml
Assay Buffer II	1 x 20ml

Relevance

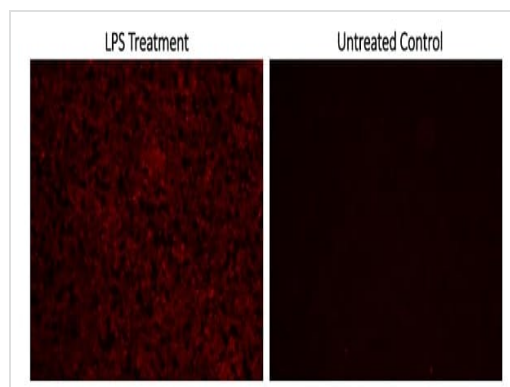
Nitric oxide (NO) is a key vertebrate biological messenger, playing an important role in neurotransmission, vascular regulation, immune responses and apoptosis. NO, also known as "endothelium-derived relaxing factor" or "EDRF", is synthesized from L-arginine, oxygen and NADPH by various NO synthases. Most of the NO in the cell is oxidized to nitrite (NO_2^-) and nitrate (NO_3^-), and therefore the concentrations of these anions are generally as a quantitative measure of NO production.

Images



Nitric Oxide Assay Kit (Fluorometric-Orange)
(ab219932)

Exogenous nitric oxide (NO) production upon DEA/NONOate treatment (NO donor). CHO-K1 (blue) and HeLa (red) cells were incubated with NO stain working solution at 37°C for 30 minutes, and then removed to stop the staining. Cells were further treated with or without 1mM DEA/NONOate in HBSS buffer (with 10 mM HEPES (pH=6.2)) at 37°C for 30 minutes. The solution in each well was removed and Assay Buffer II was added before measuring fluorescence. The fluorescence signal was monitored at Ex/Em = 540/590 nm (cut off = 570 nm) with bottom read mode using a FlexStation microplate reader (Molecular Devices).



Nitric Oxide Assay Kit (Fluorometric-Orange)
(ab219932)

Endogenous nitric oxide (NO) production in RAW 264.7 macrophage cells. Raw 264.7 cells were seeded overnight (10^5 cells/well/100 μL) in a black wall/clear bottom 96-well plate. Left: cells were treated with 20 $\mu\text{g}/\text{mL}$ of lipopolysaccharide (LPS) and 1 mM L-Arginine (L-Arg) and co-incubated with 0.5X NO stain working solution at 37°C for 16 hours. Right: control RAW 264.7 cells were left untreated in cell culture medium and co-incubated with 0.5X NO stain working solution at 37°C for 16 hours. The solution in each well was removed and Assay Buffer II was added before fluorescence measurement. The fluorescence signal was measured using fluorescence microscope with a TRITC filter.

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.com/abpromise> or contact our technical team.

Terms and conditions

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