

Peroxynitrite Assay Kit (Cell-based, Flow cytometry)
ab233470

1 References 1 Image

Overview

Product name	Peroxynitrite Assay Kit (Cell-based, Flow cytometry)
Detection method	Fluorescent
Assay type	Cell-based (quantitative)
Product overview	Due to its extremely short half-life and low steady-state concentration, it has been challenging to detect and understand the role of peroxynitrite (ONOO ⁻) in biological systems. In order to address this unmet need, ab233470 Peroxynitrite Assay Kit (Cell-based, Flow cytometry) provides a sensitive tool to monitor ONOO ⁻ levels in living cells. Peroxynitrite Sensor Green is developed as an excellent fluorescent probe, which can specifically react with intercellular ONOO ⁻ to generate a bright green fluorescent product. This kit is optimized for flow cytometry.

Notes	<p>Peroxynitrite (ONOO⁻) is a strong oxidizing species and a highly active nitrating agent. Peroxynitrite is formed from the reaction between superoxide radicals and nitric oxide generated in cells. It can damage a wide array of biomolecules including proteins, enzymes, lipids and nucleic acids, eventually contributing to cell death. Meanwhile, peroxynitrite can also have protective activities in vivo by contributing to host-defense responses against invading pathogens. Therefore, peroxynitrite is an essential biological oxidant involved in a broad range of physiological and pathological processes.</p>
-------	--

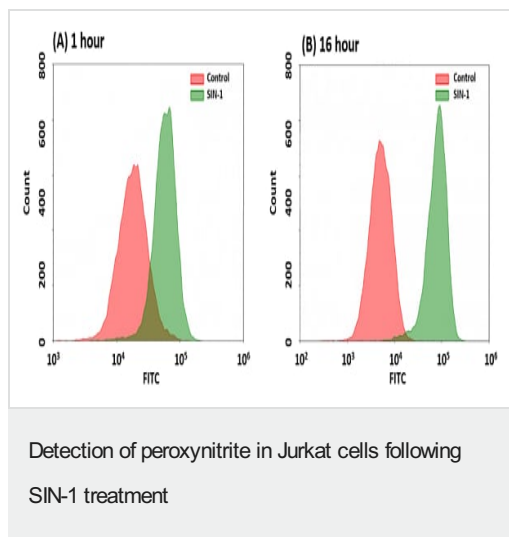
Platform	Microplate reader, Flow cytometer
----------	-----------------------------------

Properties

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

Components	100 tests
DMSO	1 x 100µl
Peroxynitrite Sensor Green	2 vials

Images



(A) Jurkat cells were co-incubated with Peroxynitrite Sensor Green and 200 μ M SIN-1 in full medium at 37 °C for 1 hour.

(B) Cells were pre-stained with Peroxynitrite Sensor Green for 1 hour, washed with PBS and then incubated with 200 μ M SIN-1 in full medium at 37 °C for 16 hours.

Cells stained with Peroxynitrite Sensor Green but without SIN-1 treatment were used as a control. Fluorescence intensity was measured using an ACEA NovoCyte flow cytometer in the FITC channel.

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