

Product datasheet

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

1 Image

Overview

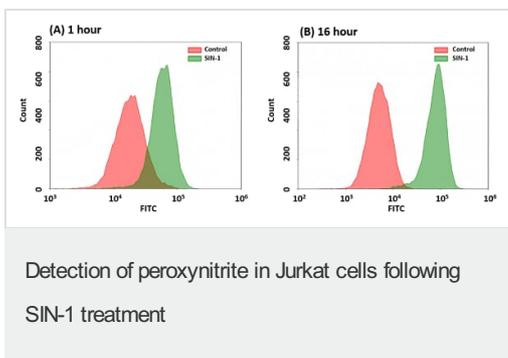
|                         |   |
|-------------------------|---|
| <b>Product name</b>     | Peroxynitrite Assay Kit (Cell-based, Flow cytometry)  |
| <b>Detection method</b> | Fluorescent   |
| <b>Assay type</b>       | Cell-based (quantitative)   |
| <b>Product overview</b> | Due to its extremely short half-life and low steady-state concentration, it has been challenging to detect and understand the role of peroxynitrite (ONOO <sup>-</sup> ) 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 <sup>-</sup> levels in living cells. Peroxynitrite Sensor Green is developed as an excellent fluorescent probe, which can specifically react with intercellular ONOO <sup>-</sup> to generate a bright green fluorescent product. This kit is optimized for flow cytometry.                |
| <b>Notes</b>            | Peroxynitrite (ONOO <sup>-</sup> ) 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. |
| <b>Platform</b>         | 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  $\mu$ M SIN-1 in full medium at 37  $^{\circ}$ C for 1 hour.

(B) Cells were pre-stained with Peroxynitrite Sensor Green for 1 hour, washed with PBS and then incubated with 200  $\mu$ M SIN-1 in full medium at 37  $^{\circ}$ 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