

Peroxynitrite Assay Kit (Fluorometric) ab233469

2 References 2 Images

Overview

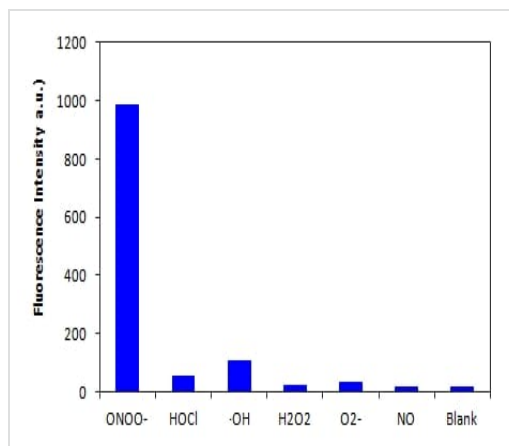
Product name	Peroxynitrite Assay Kit (Fluorometric)
Detection method	Fluorescent
Assay type	Cell-based (quantitative)
Product overview	<p>Peroxynitrite Assay Kit (Fluorometric) (ab233469) provides a sensitive tool to measure peroxynitrite in solution.</p> <p>Peroxynitrite Sensor Green reacts with <math>\text{ONOO}^-</math> to generate a bright green fluorescent product. It specifically reacts with <math>\text{ONOO}^-</math> with high selectivity over other reactive oxygen species (ROS) and reactive nitrogen species (RNS). This kit can be used with a fluorescence microplate reader and spectrometer.</p>
Notes	<p>Peroxynitrite (<math>\text{ONOO}^-</math>) is a strong oxidizing species and a highly active nitrating agent. It 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. Due to its extremely short half-life and low steady-state concentration, it has been challenging to detect and quantify peroxynitrite in solution.</p>
Platform	Microplate reader

Properties

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

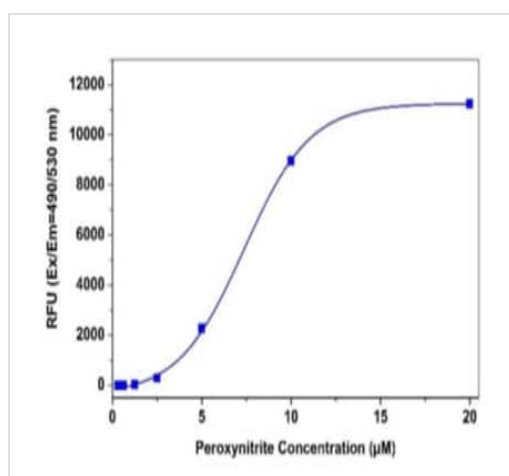
Components	100 tests
Assay Buffer	1 x 20ml
DMSO	1 x 100µl
Peroxynitrite Green reagent	1 vial

Images



Fluorescence response of Peroxynitrite Sensor  
Green to different reactive oxygen species.

The fluorescence intensities were measured with Ex/Em = 490/530 nm.



Example standard curve of Peroxynitrite showing concentration against absorbance.

Peroxynitrite was measured with the Peroxynitrite Quantification Kit (Fluorometric) (ab233469) on a solid black 96-well plate using a Gemini microplate reader. As low as 1.25 µM was detected. (Δ Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.)

**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