

Cellular ROS/RNS Assay Kit ab139473

★★★★★ 1 Abreviews 7 References

概述

产品名称	Cellular ROS/RNS Assay试剂盒
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based
种属反应性	与反应: Mammals, Other species
产品概述	Cellular ROS/RNS Assay Kit (ab139473) is designed to directly monitor real time reactive oxygen and/or nitrogen species (ROS/RNS) production in live cells using fluorescence microscopy. The kit includes three fluorescent dye reagents: NO Detection Reagent (Red), Oxidative Stress Detection Reagent (Green) for total ROS detection, and Superoxide Detection Reagent (Orange). It also includes positive controls, pyocyanin and L-arginine, which are common inducers of ROS and NO production, respectively, as well as negative controls, N-acetyl-L-cysteine and c-PTIO (common scavengers of ROS and NO, respectively). Through the combination of the three detection probes with a set of specific inhibitors and activators, this kit enables discrimination among superoxide, nitric oxide and peroxynitrite.

说明	<p>Related products</p> <p>Review the oxidative stress marker and assay guide, or the full metabolism assay guide to learn about more assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also how to assay metabolic function in live cells using your plate reader.</p>
----	---

平台	Fluorescence microscope
----	-------------------------

性能

存放说明	Please refer to protocols.
------	----------------------------

组件	200 tests
10 X Wash Buffer	1 x 15ml
NO Detection Reagent (Red)	1 x 60µl
NO Inducer (L-Arginine)	1 x 100µl
NO Scavenger (c-PTIO)	1 x 400nmole

组件	200 tests
Oxidative Stress Detection Reagent (Green)	1 x 300nmole
ROS Inducer (Pyocyanin 1 μ mole)	1 vial
ROS Inhibitor (N-acetyl-L-cysteine)	2 x 10mg
Superoxide Detection Reagent (Orange)	1 x 300nmole

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.cn/abpromise> or contact our technical team.

Terms and conditions

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