

## Product datasheet

# Cellular ROS/RNS Assay Kit ab139473

★★★★☆ 1 Abreviews 2 References

### Overview

<b>Product name</b>	Cellular ROS/RNS Assay Kit
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	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.
<b>Notes</b>	<p><b>Related products</b></p> <p>Review the <a href="#">oxidative stress marker and assay guide</a>, or the full <a href="#">metabolism assay guide</a> 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>
<b>Platform</b>	Fluorescence microscope

### Properties

**Storage instructions** Please refer to protocols.

Components	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

Components	200 tests
Oxidative Stress Detection Reagent (Green)	1 x 300nmole
ROS Inducer (Pyocyanin 1 $\mu$ 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.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