

# Cellular ROS Assay Kit (Orange) ab186028

[3 References](#) [3 Images](#)

### Overview

---

|                         |   |
|-------------------------|---|
| <b>Product name</b>     | Cellular ROS Assay Kit (Orange)   |
| <b>Sample type</b>      | Adherent cells, Suspension cells  |
| <b>Product overview</b> | Cellular ROS Assay Kit (Orange) ab186028 uses an ROS sensor to quantify ROS in live cells. The orange dye used in the ROS assay protocol is cell-permeable and generates orange fluorescence when it reacts with ROS. The fluorescence signal of the dye can be measured by fluorescence microscopy, high-content imaging, microplate fluorometry, or flow cytometry. |

ab186028 provides a sensitive fluorometric, one-step assay to detect intracellular ROS (especially superoxide and hydroxyl radical) in live cells within 1 hour. The ROS assay can be performed in a convenient 96-well or 384-well microtiter-plate format using either a fluorescence microplate reader at Ex/Em = 540/570 nm or a fluorescent microscope with TRITC filter.

**Notes** Previously called Cellular Reactive Oxygen Species Detection Assay Kit (Orange Fluorescence). Reactive oxygen species (ROS) are natural byproducts of the normal metabolism of oxygen and play important roles in cell signaling. However, during oxidative stress-related states, ROS levels can increase dramatically. The accumulation of ROS results in significant damage to cell structures. The role of oxidative stress in cardiovascular disease, diabetes, osteoporosis, stroke, inflammatory diseases, a number of neurodegenerative diseases and cancer has been well established. The ROS measurement will help to determine how oxidative stress modulates varied intracellular pathways.

#### Related products

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.

To measure reactive oxygen species within cells, we recommend [DCFDA / H2DCFDA - Cellular ROS Assay Kit ab113851](#). Alternative ROS assays are available in orange (ab186028), red ([ab186027](#)), and deep red ([ab186029](#)). [ab238535](#) is used to measure ROS in biofluids, culture supernatants and cell lysates.

For assays designed to differentiate ROS, superoxides, and reactive nitrogen species: to assay ROS and superoxides use [ab139476](#); to assay ROS, superoxides, and reactive nitrogen species use [ab139473](#); to assay superoxides use [ab219943](#).

**Platform** Microplate reader, Fluor. microscope, Flow cyt.

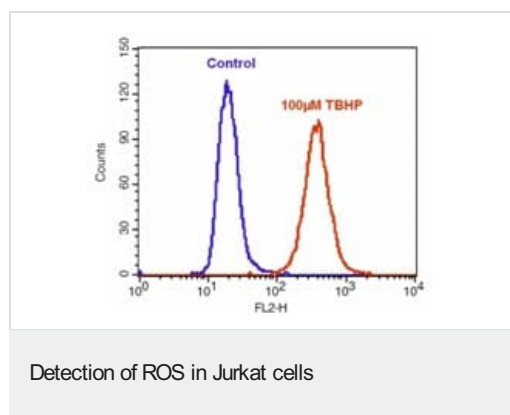
## Properties

### Storage instructions

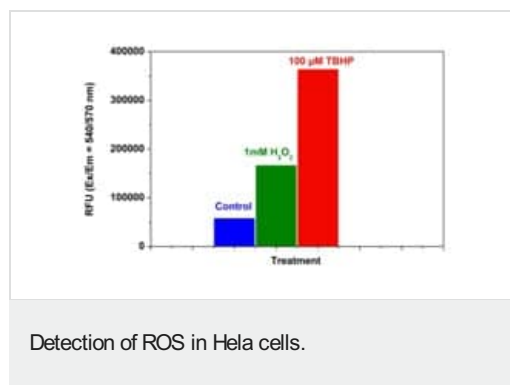
Store at -20°C. Please refer to protocols.

| Components                   | 200 tests |
|------------------------------|-----------|
| Assay Buffer                 | 1 x 20ml  |
| DMSO                         | 1 x 100µl |
| ROS Orange Dye (Lyophilized) | 1 vial    |

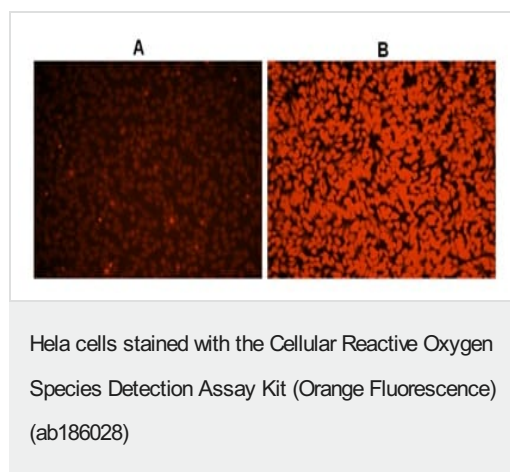
## Images



Jurkat cells were treated without (Blue) or with 100µM TBHP (Red) for 30min at 37 °C, and then loaded with ROS orange in a 5% CO<sub>2</sub>, 37 °C incubator for 1 hour. The fluorescent intensities were measured with a flow cytometer using FL2 channel



HeLa cells were seeded overnight at 15,000 cells/90 µl/well in a black wall/clear bottom 96-well plate. The cells were untreated (control) or treated with 1 mM H<sub>2</sub>O<sub>2</sub> or 100 µM TBHP for 30 minutes at 37 °C. The ROS Orange assay solution (100 µl/well) was added and incubated in a 5% CO<sub>2</sub>, 37°C incubator for 1 hour. The fluorescence signal was monitored at Ex/Em = 540/570 nm (cut off = 550 nm) with bottom read mode.



HeLa cells stained with the Cellular Reactive Oxygen Species Detection Assay Kit (Orange Fluorescence) (ab186028) in a black wall/clear bottom 96-well plate. A: Untreated control cells. B: Cells treated with 100 µM TBHP for 30min before staining.

**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