

Cell Meter™ Mitochondrial Hydroxyl Radical Detection Kit *Red Fluorescence*

Catalog number: 16055

Unit size: 200 Tests

| Component | Storage | Amount |
|-----------------------------|---|------------------|
| Component A: MitoROS™ OH580 | Freeze (<-15 °C), Minimize light exposure | 1 vial |
| Component B: Assay Buffer | Freeze (<-15 °C) | 1 bottle (50 mL) |
| Component C: DMSO | Freeze (<-15 °C) | 1 vial (100 µL) |

OVERVIEW

The detection of intracellular hydroxyl radical is of central importance to understanding proper cellular redox regulation and the impact of its dysregulation on various pathologies. The hydroxyl radical ($\cdot\text{OH}$) is one of the reactive oxygen species (ROS) highly reactive with other molecules to achieve stability. In general, hydroxyl radical is considered to be a harmful by-product of oxidative metabolism, which can cause molecular damage in living system. It shows an average lifetime of 10-9 nano seconds and can react with nearly every biomolecule such as nuclear DNA, mitochondrial DNA, proteins and membrane lipids. AAT Bioquest's Cell Meter™ Mitochondrial Hydroxyl Radical Detection Kit is optimized for detecting hydroxyl radical in mitochondria. MitoROS™ OH580 is live-cell permeant probe and can rapidly and selectively target hydroxyl radical in live cells. It generates red fluorescence when it reacts with $\cdot\text{OH}$, and can be easily read. Cell Meter™ Mitochondrial Hydroxyl Radical Detection Kit provides a sensitive fluorimetric probe to detect $\cdot\text{OH}$ in live cells with one hour incubation. This kit can be used for fluorescence microplate readers and fluorescence microscopy applications.

AT A GLANCE

Protocol summary

1. Prepare cells
2. Incubate cells with MitoROS™ OH580 working solution at 37°C for 60 minutes
3. Incubate cells with test compounds (to induce $\cdot\text{OH}$)
4. Monitor the fluorescence increase at Ex/Em = 540/590 nm

Important Thaw all the components at room temperature before use.

KEY PARAMETERS

| | |
|------------------------------|--------------------------------|
| Instrument: | Fluorescence microplate reader |
| Excitation: | 540 nm |
| Emission: | 590 nm |
| Cutoff: | 570 nm |
| Instrument specification(s): | Bottom read mode |
| Recommended plate: | Black wall/clear bottom |

| | |
|--------------------|-------------------------|
| Instrument: | Fluorescence microscope |
| Excitation: | Cy3/TRITC filter set |
| Emission: | Cy3/TRITC filter set |
| Recommended plate: | Black wall/clear bottom |

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles.

1. MitoROS™ OH580 stock solution (500X):

Add 50 µL of DMSO (Component C) into the vial of MitoROS™ OH580 (Component A), and mix them well.

Note 25 µL of stock solution is enough for 1 plate.

Note Unused portion can be aliquoted and stored at $\leq -20^{\circ}\text{C}$ for more than one month if the tubes are sealed tightly and kept from light. Avoid repeated freeze-

thaw cycles.

PREPARATION OF WORKING SOLUTION

Add 25 µL of 500X DMSO reconstituted MitoROS™ OH580 stock solution into 10 mL of Assay Buffer (Component B). Mix well.

Note This working solution is stable for at least 2 hours at room temperature.

PREPARATION OF CELL SAMPLES

For guidelines on cell sample preparation, please visit <https://www.aatbio.com/resources/guides/cell-sample-preparation.html>

SAMPLE EXPERIMENTAL PROTOCOL

1. Remove medium, and add 100 µL/well (96-well plate) or 25 µL/well (384-well plate) of MitoROS™ OH580 working solution into the cell plate. Incubate cells at 37°C for 60 minutes.
2. To induce hydroxyl radical, treat cells with test compounds in your desired buffer (such as PBS or HBBS) at 37°C for a desired period of time, protected from light.

Note We treated HeLa cells with Fenton reaction (10 µM CuCl_2 and 100 µM H_2O_2) at 37°C for 1 hour to induce exogenous hydroxyl radical. See Figure 1 for details. We treated RAW 264.7 cells with PMA (phorbol 12-myristate 13-acetate) in growth medium at 37°C for 4 hours to stimulate endogenous hydroxyl radical.
3. Wash cells 2 - 3 times with HBBS or DPBS, and add 100 µL Assay Buffer (Component B) to each well.
4. Monitor the fluorescence signal in cells using fluorescence microscope with a TRITC filter set, or measure fluorescence increase using fluorescence microplate reader at Ex/Em = 540/590 nm (cut off = 570 nm) with bottom read mode.

EXAMPLE DATA ANALYSIS AND FIGURES

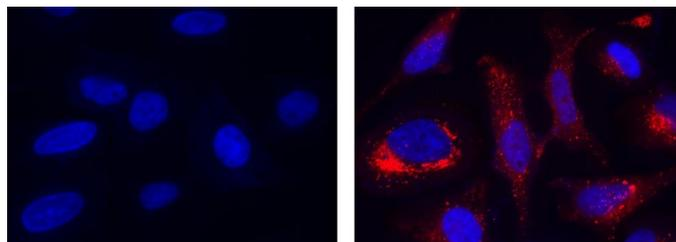


Figure 1. Fluorescence images of hydroxyl radical measurement in HeLa cells using MitoROS™ OH580 (Cat#16055). HeLa cells were incubated with MitoROS™ OH580 working solution at 37 °C for 1 hour, then washed once with HBBS. Fenton Reaction: Cells were then treated with 10 µM CuCl_2 and 100 µM H_2O_2 in 1X HBBS buffer at 37 °C for 1 hour. Control: HeLa cells were kept in 1X HBBS buffer without

treatment. After washing 3 times with HHBS, HeLa cells were measured using a fluorescence microscope with a TRITC filter set (Red). Cell nuclei were stained with Hoechst 33342 (Cat#17530, Blue).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.