

Cell Meter™ Fluorimetric Intracellular Peroxynitrite Assay Kit

Green Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 16315 (100 assays)	Keep in freezer Protect from light	Fluorescence microplate readers, Fluorescence Microscope

Introduction

Peroxynitrite (ONOO⁻) is a strong oxidizing species and a highly active nitrating agent. Peroxynitrite is formed from the reaction between superoxide radicals and nitric oxide generated in cells. It can cause damages to a wide array of biomolecules including proteins, enzymes, lipids and nucleic acids, eventually contributing to cell death. Meanwhile, peroxynitrite can also have protective activities *in vivo* by contributing to host-defense responses against invading pathogens. Therefore, peroxynitrite is an essential biological oxidant involved in a board range of physiological and pathological processes. Due to its extremely short half-life and low steady-state concentration, it has been challenging to detect and understand the role of peroxynitrite in biological systems. AAT Bioquest's Cell Meter™ Fluorimetric Intracellular Peroxynitrite (ONOO⁻) Assay Kit has been developed to address this unmet need. It provides a sensitive tool to monitor ONOO⁻ level in living cells. AAT Bioquest's DAX-J2™ PON Green is developed as an excellent fluorescent probe, which can specifically react with intercellular ONOO⁻ to generate a bright green fluorescent product. This kit can be used in fluorescence imaging and fluorescence microplate reader.

Kit Components

Components	Amount
Component A: DAX-J2™ PON Green	1 vial
Component B: Assay Buffer	1 vial (1 mL/vial)
Component C: DMSO	1 vial (100 µL/vial)

Assay Protocol for One 96-Well Plate

Brief Summary

Prepare cells in growth medium → Co-incubate cells with test compounds and DAX-J2™ PON Green working solution → Monitor fluorescence intensity at Ex/Em = 490/530 nm

1. Prepare cells:

- 1.1. **For adherent cells:** Plate cells overnight in growth medium at 20,000 to 80,000 cells/well/90 µL for a 96-well plate or 5,000 to 20,000 cells/well/22.5 µL for a 384-well plate.
- 1.2. **For non-adherent cells:** Centrifuge the cells from the culture medium and suspend the cell pellets in culture medium at 80,000-200,000 cells/well/90 µL for a 96-well poly-D lysine plate or 20,000-50,000 cells/well/22.5 µL for a 384-well poly-D lysine plate. Centrifuge the plate at 800 rpm for 2 minutes with brake off prior to your experiment.
Note: Each cell line should be evaluated on an individual basis to determine the optimal cell density.

2. Prepare working solution:

- 2.1 **Prepare DAX-J2™ PON Green stock solution (500X):** Add 20 µL of DMSO (Component C) into the vial of DAX-J2™ PON Green (Component A), and mix them well.
Note: 20 µL of reconstituted DAX-J2™ PON Green stock solution is enough for 1 plate. Unused portion can be aliquoted and stored at ≤ -20 °C for more than one month if the tubes are sealed tightly and kept from light. Avoid repeated freeze-thaw cycles.
- 2.2 **Prepare DAX-J2™ PON Green working solution (10X):** Add 10 µL of 500X DMSO reconstituted DAX-J2™ Peroxynitrite Sensor stock solution (from Step 2.1) into 500 µL of Assay Buffer (Component B), and mix them well.
Note: The working solution is not stable; prepare it as needed before use.

3. Run the peroxynitrite assay:

- 3.1 Add 10 μL /well (96-well plate), or 2.5 μL /well (384-well plate) of DAX-J2™ PON Green working solution (from Step 2.2) in 90 μL (96-well plate) or 22.5 μL (384-well plate) cell culture per well in the cell plate (from Step 1).

Note: It is not necessary to wash cells before staining. It's recommended to stain the cells in full medium.

- 3.2 Co-incubate cells with DAX-J2™ PON Green with test compounds in full medium or in your desired buffer at 37°C for desired period of time, protected from light. For control wells (untreated cells), add the corresponding amount of compound buffer.

Note 1: It's recommended to stain the cells in full medium. However, if tested compounds are serum sensitive, growth medium and serum factors can be aspirated away before staining. Add 90 μL /well (96-well plate) and 22.5 μL /well (384-well plate) of 1X Hank's salt solution and 20 mM Hepes buffer (HHBS) or the buffer of your choice after aspiration. Alternatively, cells can be stained in serum-free media.

Note 2: We co-incubated RAW 264.7 macrophage cells with 50-200 μM SIN-1 and DAX-J2™ PON Green in full medium at 37°C for 1 hour to induce peroxynitrite. See Figure 1 for details.

- 3.3 Alternatively, stain cells with DAX-J2™ PON Green at 37°C for 1 hour, protected from light (as in Step 3.1). Remove the working solution, then treat cells with test compounds in full medium or in your desired buffer at 37°C for desired period of time.
- 3.4 Monitor the fluorescence increase using microplate reader at Ex/Em = 490/530 nm (cut off = 515 nm) with bottom read mode, or take images using fluorescence microscope with a FITC filter.

Data Analysis

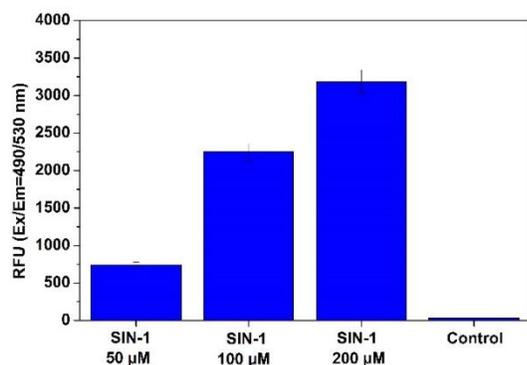


Figure 1. Detection of peroxynitrite in living cells upon SIN-1 treatment using Cell Meter™ Fluorimetric Intracellular Peroxynitrite Assay Kit (Cat#16315). RAW 264.7 cells at 100,000 cells/well/100 μL were seeded overnight in a Costar black wall/clear bottom 96-well plate. Cells were co-incubated with DAX-J2™ PON Green working solution and SIN-1 at the concentration from 50 to 200 μM at 37 °C for 1 hour. Cells incubated with DAX-J2™ PON Green without SIN-1 treatment were used as control. The fluorescence signal were monitored at Ex/Em = 490/530 nm (cut off = 515 nm) with bottom read mode using a FlexStation microplate reader (Molecular Devices).

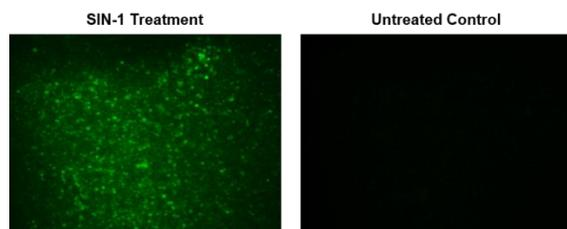


Figure 2. Fluorescence images of intracellular peroxynitrite in RAW 264.7 macrophage cells using Cell Meter™ Fluorimetric Intracellular Peroxynitrite Assay Kit (Cat#16315). Raw 264.7 cells at 100,000 cells/well/100 μL were seeded overnight in a Costar black wall/clear bottom 96-well plate. SIN-1 Treatment: Cells were co-incubated with DAX-J2™ PON Green and 100 μM SIN-1 at 37 °C for 1 hour. Untreated control: The RAW 264.7 cells were incubated with DAX-J2™ PON Green without SIN-1 treatment. The fluorescence signals were measured using a fluorescence microscope with a FITC filter.

References

1. Beckman JS and Koppenol WH. 1996. Nitric oxide, superoxide, and peroxynitrite: the good, the bad, and ugly. *Am J Physiol.* 271:1424-1437.
2. Radi R. 2013. Peroxynitrite, a stealthy biological oxidant. *JBC Papers in Press.* 288:26464-26472.
3. Szabó C, Ischiropoulos H and Radi R. 2007. Peroxynitrite: biochemistry, pathophysiology and development of therapeutics. *Nature Reviews Drug Discovery.* 6:662-680.
4. Prolo C, Alvarez MN and Radi R. 2014. Peroxynitrite, a potent macrophage-derived oxidizing cytotoxin to combat invading pathogens. *Biofactors.* 40:215-225.