

# Amplite™ Fluorimetric Peroxynitrite Quantification Kit \*Green Fluorescence\*

Catalog number: 16316  
Unit size: 100 Tests

Component	Storage	Amount
Component A: DAX-J2™ PON Green 99	Freeze (<-15 °C), Dessicated, Avoid Light	1 vial
Component B: Assay Buffer	Freeze (<-15 °C), Avoid Light	1 bottle (20 mL)
Component C: DMSO	Freeze (<-15 °C), Avoid Light	1 vial (100 µL)

## OVERVIEW

Peroxynitrite (ONOO<sup>-</sup>) 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 damage a wide array of biomolecules including proteins, enzymes, lipids and nucleic acids, eventually contributing to cell death. Due to its extremely short half-life and low steady-state concentration, it has been challenging to detect and quantify peroxynitrite in solution. In order to address this need, AAT Bioquest's Amplite™ Fluorimetric Peroxynitrite Quantification Kit provides a sensitive tool to measure ONOO<sup>-</sup> in solution. DAX-J2™ PON Green 99 reacts with ONOO<sup>-</sup> to generate a bright green fluorescent product. It specifically reacts with ONOO<sup>-</sup> with high selectivity over other reactive oxygen species (ROS) and reactive nitrogen species (RNS). This kit can be used with a fluorescence microplate reader and spectrometer.

## AT A GLANCE

**Important** Thaw all the kit components to room temperature before use.

## KEY PARAMETERS

Instrument:	Fluorescence microplate reader
Excitation:	490 nm
Emission:	530 nm
Cutoff:	515 nm
Recommended plate:	Solid black

## 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. DAX-J2™ PON Green 99 stock solution (500X):

Add 20 µL of DMSO (Component C) into the vial of DAX-J2™ PON Green 99 (Component A) to make 500X stock solution.

**Note** 20 µL of reconstituted DAX-J2™ PON Green 99 stock solution is enough for 1 plate. Keep from light.

### 2. Peroxynitrite (ONOO<sup>-</sup>) stock solution (not provided):

Peroxynitrite stock solution was synthesized according to literature report. Briefly, a mixture of sodium nitrite (0.6 M) and hydrogen peroxide (0.7 M) was acidified with hydrochloric acid (0.6 M), and sodium hydroxide (1.5 M) was added within 1 - 2 seconds to make the solution alkaline. The excess hydrogen peroxide was removed by passing the solution through a short column of manganese dioxide. The extinction coefficient of peroxynitrite solution in 0.1 M NaOH is 1670 M<sup>-1</sup>cm<sup>-1</sup> at 302 nm. The ONOO<sup>-</sup> stock solution is not stable; we highly recommend make it fresh to use.

## PREPARATION OF STANDARD SOLUTION

### Peroxynitrite standard

For convenience, use the Serial Dilution Planner:

<https://www.aatbio.com/tools/serial-dilution/16316>

Dilute Peroxynitrite (ONOO<sup>-</sup>) stock solution in Assay buffer (Component B) to have 20 µM ONOO<sup>-</sup> standard solution, and then perform 1:2 serial dilutions to get serially diluted ONOO<sup>-</sup> standard solution (O7 - O1).

## PREPARATION OF WORKING SOLUTION

Add 20 µL of 500X DAX-J2™ PON Green 99 stock solution into 10 mL of Assay Buffer (Component B) and mix well.

**Note** This assay mixture is enough for one 96-well plate. Protect from light.

## SAMPLE EXPERIMENTAL PROTOCOL

**Table 1.** Layout of ONOO<sup>-</sup> standards and test samples in a solid black 96-well microplate. O = ONOO<sup>-</sup> Standards (O1 - O7, 0.313 to 20 µM); BL = Blank Control; TS = Test Samples.

BL	BL	TS	TS
O1	O1	...	...
O2	O2	...	...
O3	O3		
O4	O4		
O5	O5		
O6	O6		
O7	O7		

**Table 2.** Reagent composition for each well.

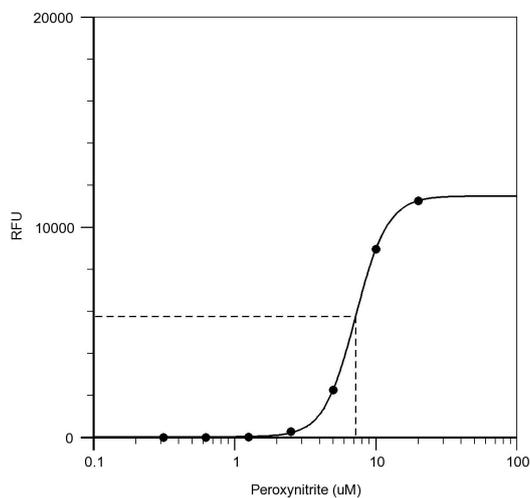
Well	Volume	Reagent
O1 - O7	50 µL	Serial Dilutions (0.313 to 20 µM)
BL	50 µL	assay buffer
TS	50 µL	test sample

1. Prepare ONOO<sup>-</sup> standards (O), blank controls (BL), and test samples (TS) according to the layout provided in Tables 1 and 2. For a 384-well plate, use 25 µL of reagent per well instead of 50 µL.
2. Add 50 µL of working solution to each well of ONOO<sup>-</sup> standard, blank control, and test samples to make the total ONOO<sup>-</sup> assay volume of 100 µL/well. For a 384-well plate, add 25 µL of working solution into each well instead, for a total volume of 50 µL/well.
3. Incubate the reaction at room temperature for 5 to 10 minutes, protected from light.
4. Monitor the fluorescence increase at Ex/Em = 490/530 nm (cutoff at 515 nm) with a fluorescence plate reader.

**EXAMPLE DATA ANALYSIS AND FIGURES**

The reading (RFU) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the baseline corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate Peroxynitrite samples. We recommend using the Online Four Parameter Logistics Calculator which can be found at:

<https://www.aatbio.com/tools/four-parameter-logistic-4pl-curve-regression-online-calculator/>



**Figure 1.** Peroxynitrite was measured with Amplite™ Fluorimetric Peroxynitrite Quantification Kit on a solid black 96-well plate using a Gemini microplate reader (Molecular Devices).

**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.