

## Amplite™ Fluorimetric Hydrogen Peroxide Assay Kit

### \*Red Fluorescence\*

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 11501 (500 assays)	Keep in freezer Avoid exposure to light	Fluorescence microplate readers

### Introduction

Hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) is a reactive oxygen metabolic by-product that serves as a key regulator for a number of oxidative stress-related states. It is involved in a number of biological events that have been linked to asthma, atherosclerosis, diabetic vasculopathy, osteoporosis, a number of neurodegenerative diseases and Down's syndrome. Perhaps the most intriguing aspect of H<sub>2</sub>O<sub>2</sub> biology is the recent report that antibodies have the capacity to convert molecular oxygen into hydrogen peroxide to contribute to the normal recognition and destruction processes of the immune system. Measurement of this reactive species will help to determine how oxidative stress modulates a variety of intracellular pathways.

This Amplite™ Fluorimetric Hydrogen Peroxide Assay Kit uses our non-fluorescent Amplite™ Red peroxidase substrate to quantify hydrogen peroxide in solutions and cell extracts. It can also be used to detect a variety of oxidase activities through enzyme-coupled reactions. The kit is an optimized “mix and read” assay that is compatible with HTS liquid handling instruments. It provides a sensitive, one-step fluorometric assay to detect as little as 3 picomoles of H<sub>2</sub>O<sub>2</sub> in a 100 µL assay volume (30 nM, Figure 1). The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and readily adapted to automation. Its signal can be easily read by either a fluorescence microplate reader at Ex/Em = ~540/590 nm or an absorbance microplate reader at ~570 nm.

#### Kit Key Features

<b>Broad Application:</b>	Can be used for quantifying hydrogen peroxide in solutions, in cell extracts and in live cells; and can also be used for detecting a variety of oxidase activities through enzyme-coupled reactions.
<b>Sensitive:</b>	Detect as low as 10 picomoles of H <sub>2</sub> O <sub>2</sub> in solution.
<b>Continuous:</b>	Easily adapted to automation without a separation step.
<b>Convenient:</b>	Formulated to have minimal hands-on time. No wash is required.
<b>Non-Radioactive:</b>	No special requirements for waste treatment.

### Kit Components

Components	Amount
Component A: Amplite™ Red Peroxidase Substrate	1 vial
Component B: H <sub>2</sub> O <sub>2</sub>	1 vial (3% stabilized solution, 200 µL)
Component C: Assay Buffer	1 bottle (100 mL)
Component D: Horseradish Peroxidase	1 vial (20 units)
Component E: DMSO	1 vial (1 mL)

## Assay Protocol for One 96-well Plate

### Brief Summary

**Prepare H<sub>2</sub>O<sub>2</sub> reaction mixture (50 µL) → Add H<sub>2</sub>O<sub>2</sub> standards or test samples (50 µL) → Incubate at room temperature for 10-30 minutes → Monitor fluorescence intensity at Ex/Em = 540/590 nm**

*Note: Thaw all the kit components at room temperature before starting the experiment.*

#### 1. Prepare stock solutions:

1.1 100X Amplite™ Red peroxidase substrate stock solution: Add 250 µL of DMSO (Component E) into the vial of Amplite™ Red Substrate (Component A). The stock solution should be used promptly; any remaining solution should be aliquoted and refrozen at -20 °C.

*Note: Avoid repeated freeze-thaw cycles and protect from light.*

1.2 20 U/mL Peroxidase stock solution: Add 1 mL of Assay Buffer (Component C) into the vial of Horseradish Peroxidase (Component D).

*Note: The unused HRP solution should be divided into single use aliquots and stored at -20 °C.*

1.3 20 mM H<sub>2</sub>O<sub>2</sub> stock solution: Add 22.7 µL of 3% H<sub>2</sub>O<sub>2</sub> (0.88 M, Component B) into 977 µL of Assay Buffer (Component C).

*Note: The diluted H<sub>2</sub>O<sub>2</sub> solution is not stable. The unused portion should be discarded.*

#### 2. Prepare H<sub>2</sub>O<sub>2</sub> reaction mixture:

Prepare the H<sub>2</sub>O<sub>2</sub> reaction mixture according to the following table and keep from light:

**Table 1** H<sub>2</sub>O<sub>2</sub> Reaction mixture for one 96-well plate (2X)

Components	Volume
Amplite™ Red Peroxidase Substrate Stock Solution (100X, from Step 1.1)	50 µL
20 U/ml Peroxidase Stock Solution (from Step 1.2)	200 µL
Assay Buffer (Component C)	4.75 mL
Total volume	5 mL

#### 3. Prepare serial dilutions of H<sub>2</sub>O<sub>2</sub> standard (0 to 10 µM):

**Warning 1:** The component A is unstable in the presence of thiols such as DTT and β-mercaptoethanol. Thiols higher than 10 µM (final concentration) would significantly decrease the assay dynamic range.

**Warning 2:** NADH and glutathione (reduced form: GSH) may interfere with the assay.

3.1 Add 1 µL of 20 mM H<sub>2</sub>O<sub>2</sub> solution (from Step 1.3) into 1999 µL of Assay Buffer (Component C) to get a 10 µM H<sub>2</sub>O<sub>2</sub> standard.

3.2 Take 200 µL of 10 µM H<sub>2</sub>O<sub>2</sub> standard to perform 1:3 serial dilutions to get 3, 1, 0.3, 0.1, 0.03, 0.01 and 0 µM serial dilutions of H<sub>2</sub>O<sub>2</sub> standard.

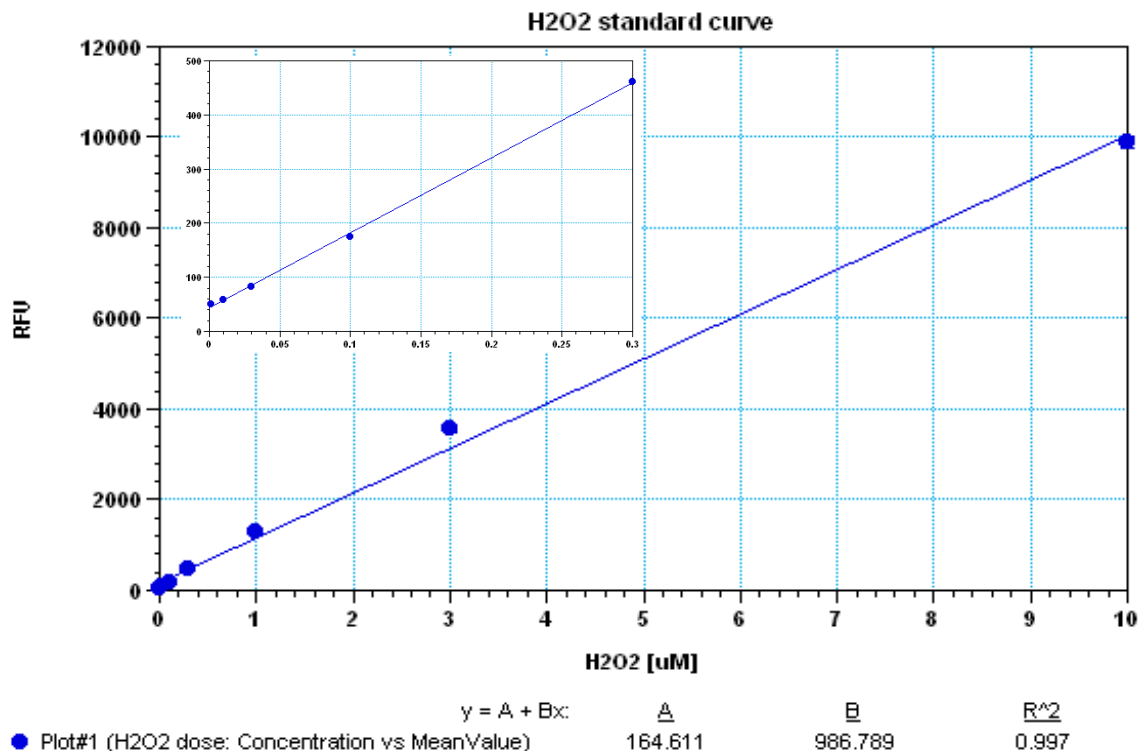
3.3 Add serial dilutions of H<sub>2</sub>O<sub>2</sub> standard and H<sub>2</sub>O<sub>2</sub>-containing test samples into a solid black 96-well microplate as described in Tables 2 and 3.



## Data Analysis

The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with the H<sub>2</sub>O<sub>2</sub> reactions. A H<sub>2</sub>O<sub>2</sub> standard curve is shown in Figure 1.

*Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.*



**Figure 1.** H<sub>2</sub>O<sub>2</sub> dose response was measured in a 384-well black plate with the Amplite™ Florimetric Hydrogen Peroxide Assay Kit using a Gemini fluorescence microplate reader (Molecular Devices). As low as 0.03 µM H<sub>2</sub>O<sub>2</sub> can be detected with 30 minutes incubation (n=3). The insert shows the low levels of H<sub>2</sub>O<sub>2</sub> detection.

## References

- Hoffmann, O. M., Becker, D., and Weber, J. R. (2007) *J Cereb Blood Flow Metab.*
- Funk, R. S., and Krise, J. P. (2007) *Mol Pharm* 4, 154-9.
- Krebs, B., Wiebelitz, A., Balitzki-Korte, B., Vassallo, N., Paluch, S., Mitteregger, G., Onodera, T., Kretschmar, H. A., and Herms, J. (2007) *J Neurochem* 100, 358-67.
- Yang, Y., Xu, S., An, L., and Chen, N. (2007) *J Plant Physiol.*
- Lee, J. E., Kim, H., Jang, H., Cho, E. J., and Youn, H. D. (2007) *J Neurochem.*
- Goth, L. (2006) *Redox Rep* 11, 281-2.
- Laloi, C., Stachowiak, M., Pers-Kamczyc, E., Warzych, E., Murgia, I., and Apel, K. (2007) *Proc Natl Acad Sci U S A* 104, 672-7.

**Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.**