



SensoLyte[®] ADHP Hydrogen Peroxide Assay Kit **Fluorimetric**

Revision# 1.2	Last Updated: July 2021
Catalog #	AS-71112
Kit Size	500 Assays (96-well plate)

- **Convenient Format:** Complete kit includes all the assay components.
- **Optimized Performance:** Optimal conditions for quantifying hydrogen peroxide and detecting oxidase.
- **Enhanced Value:** Less expensive than the sum of individual components.
- **High Speed:** Minimal hands-on time.
- **Assured Reliability:** Detailed protocol and references are provided.

Kit Components, Storage and Handling

Component	Description	Quantity
Component A	ADHP	10 mM, 250 μ L
Component B	H ₂ O ₂ standard	1 vial
Component C	Assay buffer	60 mL
Component D	HRP, Horseradish peroxidase	5 vials, 100 μ L/vial

Other Materials Required (but not provided)

- 96-well microplate: Black, flat-bottom microplates with non-binding surface.
- Fluorescence microplate reader: Capable of detecting emission at 590 nm with excitation at 530-560 nm.

Storage and Handling

- For convenience, Component C can be stored at room temperature.

Introduction

Reactive oxygen species (ROS) play an important role in a variety of biological events, such as inflammation, ischemia and reperfusion, and neurodegeneration. Hydrogen peroxide (H_2O_2) is membrane permeable and is more stable than other ROS. It is often chosen to represent the ROS released by cell or cell organelles (e.g. mitochondria,¹ activated leukocytes²). H_2O_2 is also a co-product of many oxidase-catalyzed reactions. Consequently, it can serve as an indicator of the activity of oxidases (e.g. NADPH oxidase³, glucose oxidase⁴, and monoamine oxidase⁵).

The SensoLyte[®] ADHP Hydrogen Peroxide Assay Kit provides a convenient, highly sensitive fluorescent assay for quantifying H_2O_2 in solutions, in cell extracts and in live cells. In the enzyme-coupled reaction, non-fluorescent ADHP (10-Acetyl-3, 7-dihydroxyphenoxazine) can be oxidized to the strongly fluorescent resorufin in presence of H_2O_2 and horseradish peroxidase (HRP). The signal of resorufin can be easily read by a fluorescence microplate reader at Ex/Em=530-560 nm/590 nm.

Protocol

Note: Warm all kit components to room temperature before starting the experiment.

1. Prepare stock solution.

- H_2O_2 stock solution (1 M): Add 100 μL of deionized water into the H_2O_2 vial (Component B) to get 1 M stock solution. Store this stock solution tightly capped at 4°C.

2. Set up the H_2O_2 standard curve (Optional).

- Dilute 1 M H_2O_2 stock solution to 40 μM in assay buffer (Component C). Perform 2-fold serial dilutions with the assay buffer to get 20, 10, 5, 2.5, 1.25, and 0.63 μM H_2O_2 solutions. Add 50 μL /well of the serially diluted H_2O_2 solution. Include a negative control that does not contain any H_2O_2 .

3. Prepare test samples.

- Add 50 μL /well of samples (e.g. mitochondria¹, activated leukocytes², monoamine oxidase with its substrate benzylamine³).

Note: Extremely large amount of H_2O_2 (e.g. >100 μM) may further convert fluorescent resorufin to non-fluorescent resazurin and lead to reduction of fluorescence signal. It is necessary to test your sample with several different dilutions.

4. Prepare ADHP reaction mixture.

- Prepare fresh ADHP reaction mixture according to the following Table 1 and keep away from light.

Table 1. ADHP reaction mixture for one 96-well plate (100 assays)

Components	Volume
ADHP (Component A)	50 μL
HRP (Component D)	100 μL
Assay buffer (Component C)	4.85 mL
Total volume	5 mL

Note 1: This reaction mixture can detect 0.1 nmol of H₂O₂ with a linear range of up to 2 nmol (Figure 1). Lowering the ADHP concentration in the reaction mixture can decrease background and increase assay sensitivity. 10 μM ADHP can detect as low as 2 pmol of H₂O₂.² 2 μM ADHP was used to detect H₂O₂ produced by mitochondria¹.

Note 2: You may change the assay buffer to any buffer appropriate for your samples. For example, you may use Krebs-Ringer phosphate for detecting H₂O₂ released from activated human leukocytes² or modified buffer for mitochondria¹. You may also add stimulating reagents in the reaction mixture.²

5. Detect H₂O₂.

5.1 Add 50 μL/well of ADHP reaction mixture. Mix the reagents by gently shaking the plate for 30 sec.

5.2 Incubate the reaction at the desired temperature for 15-30 min. Measure emission at 590 nm with excitation at 530-560 nm.

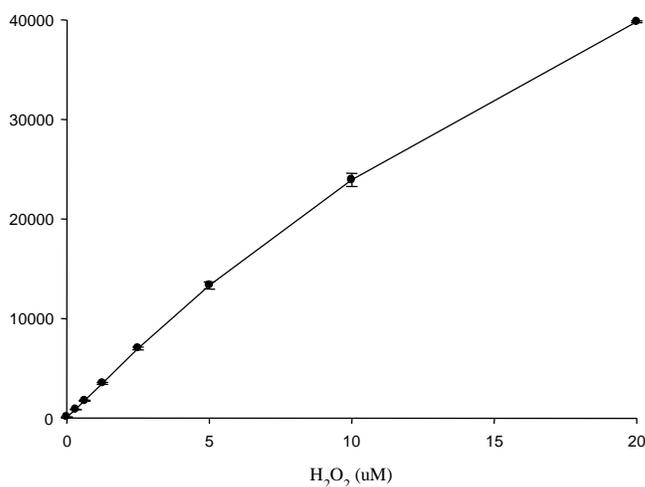


Figure 1. The standard curve of H₂O₂. H₂O₂ was serially diluted and detected according to the above protocol. With the total assay volume of 100 μL, the assay can detect as low as 1 μM (0.1 nmol) H₂O₂ with a linear range up to 20 μM (2 nmol) ($R^2 > 0.98$). (n=2, mean±S.D.)

References

1. Votyakova, T.V. and Reynolds I.J., *J. Neurochem.* **79**, 266 (2001)
2. Mohanty, J.G. et al. *J. Immunol. Methods.* **202**, 133 (1997)
3. Zhou, M. et al. *Anal. Biochem.* **253**, 162 (1997)
4. Sanchez, F.A. et al. *Anal. Biochem.* **187**, 129 (1990)
5. Youdim, M.B. and Tenne M., *Methods. Enzymol.* **142**, 617 (1987)