

Amplite™ Fluorimetric Catalase Assay Kit

Red Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 11306 (200 assays)	Keep in freezer. Avoid exposure to light.	Fluorescence microplate readers

Introduction

Catalase is a common antioxidant heme-containing redox enzyme found in nearly all living organisms that are exposed to oxygen. The enzyme is concentrated in the peroxisome subcellular organelles. Hydrogen peroxide is an ROS that is a toxic product of normal aerobic metabolism and pathogenic ROS production involving oxidase and superoxide dismutase reactions. By preventing the excessive buildup of H₂O₂, catalase allows important cellular processes which produce H₂O₂ as a by-product to take place safely.

The Amplite™ Fluorimetric Catalase Assay Kit provides a quick and sensitive method for the measurement of catalase activity. It can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. Catalase reacts with H₂O₂ to produce water and oxygen (O₂). Amplite™ Red also reacts with H₂O₂ to generate a red fluorescent product. Therefore the reduction in fluorescence intensity is proportional to catalase activity. The Amplite™ Red substrate used in the assay enables a dual recordable mode. The fluorescent signal can be easily read by either a fluorescence microplate reader at Ex/Em = 540/590 nm or an absorbance microplate reader at ~576 nm. With the Amplite™ Fluorimetric Catalase Assay Kit, we have detected as little as 30 mU/mL catalase in a 100 µL reaction volume.

Kit Key Features

Sensitive:	Detect as low as 30 mU/mL catalase.
Continuous:	Easily adapted to automation without a separation step.
Convenient:	Formulated to have minimal hands-on time.
Non-Radioactive:	No special requirements for waste treatment.

Kit Components

Components	Amount
Component A: Amplite™ Red	1 vial
Component B: H ₂ O ₂	1 vial (3% stabilized solution, 200 µL)
Component C: Assay Buffer	1 bottle (20 mL)
Component D: Horseradish Peroxidase	1 vial (20 units)
Component E: Catalase Standard	1 vial (1000 U/mL, 50uL)
Component F: DMSO	1 vial (200 µL)

Assay Protocol for One 96-well Plate

Brief Summary

Prepare catalase standards and/or test samples (50 µL) → Add H₂O₂ Assay Buffer (50 µL) → Incubate at room temperature for 10-30 minutes → Add Catalase Assay Mixture (50 µL) → Incubate at room temperature for 10-30 minutes → Monitor fluorescence increase at Ex/Em = 540/590 nm

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

Table 3. Assay mixture for one 96-well plate

Components	Volume
Amplite™ Red substrate stock solution (200X, from Step 1.1)	25 μ L
HRP stock solution (100U/mL, from Step 1.2)	15 μ L
Assay Buffer (Component C)	5.0 mL

4.4 Add 50 μ L of assay mixture (from Step 4.3) into each well of catalase standard, blank control, and test samples (see Step 3.3) to make the total assay volume of 150 μ L/well.

Note: For a 384-well plate, add 25 μ L of assay mixture into each well.

4.5 Incubate the reaction at room temperature for 15 to 30 minutes, protected from light.

4.6 Monitor the fluorescence increase at Ex/Em = 540 \pm 10/590 \pm 10 nm (optimal Ex/Em = 540/590) using a fluorescence plate reader.

Note: The contents of the plate can also be transferred into a white clear bottom plate and read by an absorbance microplate reader at the wavelength of 576 \pm 5 nm. The absorption detection has lower sensitivity compared to fluorescence reading.

Data analysis

The fluorescence intensity of blank wells (no catalase, with the assay buffer only) is used as a control. For each catalase dose, the fluorescence intensity is reported as the **difference** of the observed fluorescence intensity subtracted from that of a no-catalase control. The catalase standard curve is shown in Figure 1.

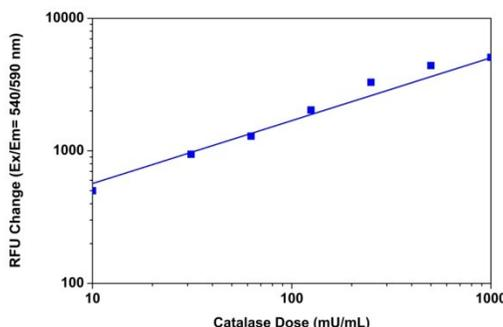


Figure 1 Catalase dose response was measured with Amplite™ Fluorimetric Catalase Assay Kit in a solid black 96-well plate using a Gemini fluorescence microplate reader (Molecular Devices). As low as 30 mU/mL catalase can be detected with 30 minutes incubation (n=3).

References

1. Porstmann, B., Porstmann, T., Nugel, E. and Evers, U. (1985). Which of the commonly used marker enzymes gives the best results in colorimetric and fluorimetric enzyme immunoassays: horseradish peroxidase, alkaline phosphatase, β -galactosidase? *J. Immunol. Meth.* **79**, 27-37.
2. Wordinger, R.J., Miller, G.W. and Nicodemus, D.S. (1987). *Manual of Immunoperoxidase Techniques, 2nd Edition*. Chicago: American Society of Clinical Pathologists Press, pp. 23-24.
3. Yolken, R.H. (1982). Enzyme immunoassays for the detection of infectious antigens in body fluids: current limitations and future prospects. *Rev. Infect. Dis.* **4(1)**, 35-68.

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 the 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.