



Hydrogen Peroxide Cell-Based Assay Kit

Item No. 600050

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

GENERAL INFORMATION	3	Materials Supplied
	4	Safety Data
	4	Precautions
	5	If You Have Problems
	5	Storage and Stability
	5	Materials Needed but Not Supplied
INTRODUCTION	6	About This Assay
PRE-ASSAY PREPARATION	7	Reagent Preparation
ASSAY PROTOCOL	9	Performing the Assay
ANALYSIS	10	Calculations
	12	Performance Characteristics
RESOURCES	13	Troubleshooting
	14	References
	15	Notes
	15	Warranty and Limitation of Remedy

GENERAL INFORMATION

Materials Supplied

For best results, upon receipt of the delivered kit, remove components and store as stated below.

Item Number	Item	96 well Quantity/Size	480 well Quantity/Size	Storage
10009322	Cell-Based Assay Buffer Tablet	1 tablet	1 tablet	RT
600051	Cell-Based Assay Hydrogen Peroxide Standard	1 vial/100 μ l	1 vial/100 μ l	4°C
600052	Cell-Based Assay Hydrogen Peroxide Detector ADHP	1 vial/100 μ l	1 vial/500 μ l	-20°C
600053	Cell-Based Assay Horseradish Peroxidase	1 vial/100 μ l	1 vial/500 μ l	-20°C
600054	Cell-Based Assay Catalase	1 vial/2 mg	1 vial/10 mg	-20°C

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



WARNING: THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

Safety Data

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent *via* email to your institution.

Precautions

Please read these instructions carefully before beginning this assay.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888
Fax: 734-971-3641
Email: techserv@caymanchem.com
Hours: M-F 8:00 AM to 5:30 PM EST

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored as directed in the **Materials Supplied** section and used before the expiration date indicated on the outside of the box.

Materials Needed But Not Supplied

1. A plate reader with the capacity to measure fluorescence using an excitation wavelength between 530-560 nm and an emission wavelength of 590 nm.
2. Serum-free cell culture media appropriate for the cells used.
3. A source of pure water; glass distilled water or HPLC-grade water are acceptable.
4. A 96-well black plate.

About This Assay

Cayman's H₂O₂ Cell-Based Assay Kit provides a simple fluorometric method for the sensitive quantitation of extracellular H₂O₂ produced by cultured cells. H₂O₂ is detected using ADHP (10-acetyl-3,7-dihydroxyphenoxazine), a highly sensitive and stable probe for H₂O₂.¹ In a horseradish peroxidase catalyzed reaction, ADHP reacts with H₂O₂ with a 1:1 stoichiometry to produce highly fluorescent resorufin.² Resorufin fluorescence can be read using an excitation wavelength between 530-560 nm and an emission wavelength of 590 nm. Catalase, an H₂O₂ scavenger, is included in the kit to check specificity of the assay. Reactive oxygen species, such as hydrogen peroxide (H₂O₂) and superoxide are generated by phagocytes and participate in damaging invading microorganisms or other biologic targets.³ This kit provides a valuable method for immunologists to assess the capacity of immune cell killing using ROS.

Reagent Preparation

1. Assay Buffer Preparation

Dissolve each Cell-Based Assay Buffer Tablet (Item No. 10009322) in 100 ml of UltraPure water. This buffer should be stable for approximately one year at room temperature.

2. Stock H₂O₂ Standard/Positive Control Preparation

Due to the instability of H₂O₂, it is recommended that the concentration of the stock H₂O₂ (Item No. 600051) be assessed prior to use. To measure the stock concentration, dilute the stock solution 1:1,000 in UltraPure water and read the absorbance at 240 nm in a quartz cuvette.

The concentration (M) is equivalent to the absorbance x the dilution factor divided by 43.6. This stock can then be diluted to 1 mM to provide a 100X working stock to begin your standard curve.

To prepare the standard for use in the H₂O₂ assay: obtain eight clean test tubes and label them #1 through #8. Aliquot 990 µl of serum-free culture medium into tube #1 and 500 µl into tubes #2-#8. Transfer 10 µl of the 1 mM stock H₂O₂ into tube #1 and mix thoroughly. Serially dilute the standard by removing 500 µl from tube #1 and place into tube #2; mix thoroughly. Repeat this procedure for tube #3 to tube #7. Do not add any standard to tube #8, your blank. The final concentrations of the standard tubes are 10, 5, 2.5, 1.25, 0.625, 0.313, 0.156, and 0 µM, respectively.

If a quantitative measurement is not required, assume the standard is approximately 8.8 M and dilute it to 10 mM for the positive control. Dilute 1:1,000 immediately prior to use. This solution is not stable and should be discarded after use.

3. Catalase Solution Preparation

Reconstitute the Cell-Based Assay Catalase (Item No. 600054) in Assay Buffer as follows: 2 mg vial (96-well kit) in 500 μ l, or 10 mg vial (480-well kit) in 2.5 ml. Keep on ice and aliquot for storage at -20°C if multiple uses are anticipated. The reconstituted catalase is stable for one month at -20°C .

4. Enzyme Reaction Solution Preparation

To make 1 ml of Enzyme Reaction Solution, sufficient for use on one 96-well plate, add 100 μ l of Cell-Based Assay Hydrogen Peroxide Detector ADHP (Item No. 600052) and 100 μ l Cell-Based Assay Horseradish Peroxidase (Item No. 600053) to 800 μ l of the Assay Buffer. This Enzyme Reaction Solution is stable for up to one hour when kept on ice. Aliquot the 480-well vial into smaller aliquots and store at -20°C to avoid repeated freeze/thaw cycles.

ASSAY PROTOCOL

Performing the Assay

Procedure

1. Seed cells in a 96-well plate at a density of 10^4 - 10^5 cells/well in 100 μ l of serum-free culture medium. If desired, compounds or vehicle controls can be added to the 100 μ l of culture medium. Control wells should be included which contain medium and experimental compounds but no cells. Include extra wells for catalase controls to determine assay specificity. We recommend that each treatment be performed in triplicate.
2. Culture the cells at 37°C for 24-48 hours, or for a period of time according to your typical experimental protocol. Centrifuge the 96-well tissue culture plate at $400 \times g$ for five minutes.
3. Transfer 80 μ l of supernatant from each well of the cultured cells or from the standard tubes prepared earlier to a new 96 well black plate. Add 10 μ l of Assay Buffer (sample activity measurement) or 10 μ l of the Catalase Solution (to test assay specificity) to appropriate wells.
4. Add 10 μ l of the Enzyme Reaction Solution to each well.
5. Within five minutes, read the fluorescence intensity of each well (excitation = 530 nm; emission = 590 nm).

Calculations

Determination of the Reaction Rate

1. Calculate the average fluorescence of each standard and sample.
2. Subtract the average fluorescence of the blank (Standard tube #8) from itself and from all other standards and samples including the catalase containing samples.
3. Plot the corrected fluorescence of the standards (from step 1 above) as a function of the final H_2O_2 concentration (μM). See Figure 1 (below) for a typical standard curve.

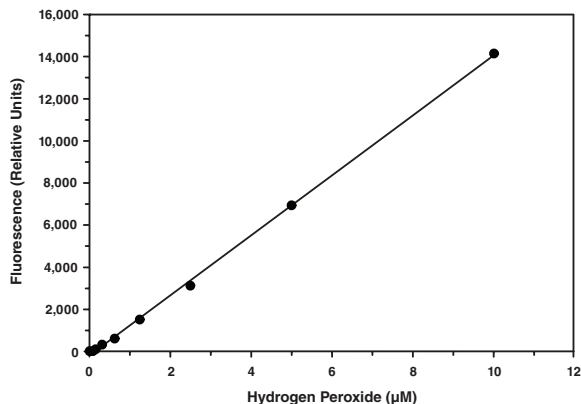


Figure 1. Example H_2O_2 standard curve

4. Subtract the catalase sample fluorescence from the non-catalase sample fluorescence to yield the corrected sample fluorescence.
5. Calculate the H_2O_2 concentration of the samples using the equation obtained from the linear regression of the standard curve substituting corrected fluorescence values for each sample.

H_2O_2 Concentration (μM) =

$$\left[\frac{\text{Corrected sample fluorescence} - (\text{y-intercept})}{\text{Slope}} \right] \times \text{Dilution}$$

If a high production of H_2O_2 in the samples is expected, serial dilution may be required to obtain values that fall on the standard curve.

Performance Characteristics

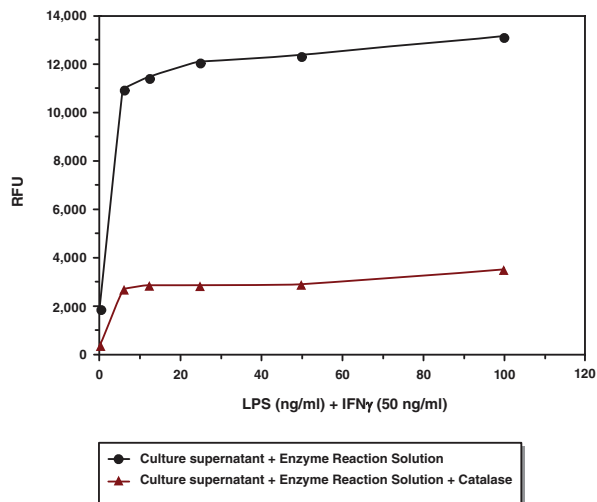


Figure 2. IFN γ and LPS together increase H₂O₂ production in RAW 264.7 cells.

RESOURCES

Troubleshooting

Problem	Possible Causes	Recommended Solutions
Erratic values; dispersion of duplicates/triplicates	A. Poor pipetting/technique B. Bubble in the well(s)	A. Be careful not to splash the contents of the wells B. Carefully tap the side of the plate with your finger to remove bubbles
No H ₂ O ₂ was detected in the sample and standard wells	Enzyme mixture was not prepared correctly	Prepare a fresh enzyme mixture and re-assay
Erratic response curve of compound treatments	Unequal number of cells in each well	Make sure each well contains the same number of cells

References

1. Zhou, M., Diwu, Z., Panchuk-Voloshina, N., *et al.* A stable nonfluorescent derivative of resorufin for the fluorometric determination of trace hydrogen peroxide: Applications in detecting the activity of phagocyte NADPH oxidase and other oxidases. *Anal. Biochem.* **253**, 162-168 (1997).
2. Amundson, D.M. and Zhou, M. Fluorometric method for the enzymatic determination of cholesterol. *J. Biochem. Biophys. Meth.* **38**, 43-52 (1999).
3. Rosen, G.M., Pou,S., Ramos, C.L. *et al.* Free radicals and phagocytic cells. *FASEB J.* **9**, 200-209 (1995).

NOTES

Warranty and Limitation of Remedy

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©07/23/2018, Cayman Chemical Company, Ann Arbor, MI, All rights reserved.
Printed in U.S.A.

