

Hydrogen peroxide colorimetric detection kit

Catalog #: ADI-907-015

96 Determination Kit

For the latest product information, including support documentation, visit us online:

www.enzolifesciences.com

Technical Support (US): 800-942-0430
Technical Support (EU): +41 61 926 8989



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research use only. Not for use in diagnostic procedures. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Description	2
Introduction	2
Precautions	2
Materials Supplied.....	3
Storage	3
Other Materials Needed	3
Sample Handling.....	4
Procedural Notes	4
Reagent Preparation	5
Assay Procedure.....	6
Calculation of Results	7
Typical Results.....	8
Typical Standard Curve.....	8
Performance Characteristics	9
Sample Recoveries	11
References.....	12
Contact Information.....	14

DESCRIPTION

The Hydrogen peroxide colorimetric detection kit is a complete kit for the quantitative determination of Hydrogen Peroxide in biological fluids and tissue culture media. Please read the complete kit insert before performing this assay. The kit is designed to measure low concentrations of H₂O₂ in biological matrices. The kit has a color reagent that contains a dye, xylene orange, in an acidic solution with sorbitol and ammonium iron sulfate that reacts to produce a purple color proportional to the concentration of H₂O₂ in the sample. The exact mechanism of the color reaction is not known, but probably involves coordinated iron reacting with H₂O₂ and the dye molecule.

INTRODUCTION

Hydrogen Peroxide (H₂O₂) is a reactive oxygen metabolic by-product that serves as a key regulator for a number of oxidative stress-related states.^{1,2} Functioning through NF kappa-B and other factors, hydroperoxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis, diabetic vasculopathy, osteoporosis, a number of neurodegenerative diseases and Down's Syndrome.³⁻¹¹ Perhaps the most intriguing aspect of H₂O₂ 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.^{12,13} Measurement of this reactive species will help to determine how oxidative stress modulates varied intracellular pathways.

PRECAUTIONS

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

1. Dispose of the contents of the plate with care. Attention should be taken in handling because of unknown effects of the contents.
2. We test this kit's performance in a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results. Refer to section titled "Sample Handling".
3. The standard is light-sensitive and should be protected from direct light for prolonged periods of time.



Handle
with care



Protect
from light

MATERIALS SUPPLIED

1. **Half-Area Microtiter Plate, 1 each, Catalog No. 80-2404:**
The plate is ready to use.
2. **Hydrogen Peroxide Standard, 0.5ml, Catalog No. 80-0941:**
A solution of Hydrogen Peroxide at 100,000ng/mL in water with preservatives.
3. **Hydrogen Peroxide Color Reagent, 11ml, Catalog No. 80-0968**
A solution of colorimetric substrate in dilute acid.
4. **Hydrogen Peroxide Assay Layout Sheet, 1 each, Catalog No. 30-0170**
5. **Plate Sealer, 2 each, Catalog No. 30-0012**



Reagents require separate storage conditions.

STORAGE

All components of this kit, except the Hydrogen Peroxide Color Reagent, are stable at 4°C until the kit's expiration date. The Hydrogen Peroxide Color Reagent must be stored at -20°C.

OTHER MATERIALS NEEDED

1. Recommended sample diluent, 50mM Phosphate, pH 6.0.
2. Precision pipets for volumes between 34µl and 1,000µl.
3. Repeater pipet for dispensing 100µl.
4. Microplate reader capable of reading between 540 and 570nm, ideally 550nm.
5. Graph paper for plotting the standard curve.

SAMPLE HANDLING

The Hydrogen peroxide colorimetric detection kit is compatible with samples in a wide range of matrices. Samples diluted sufficiently into the recommended Sample Diluent of 50mM Phosphate, pH 6.0 can be read directly from the standard curve. Please refer to the Sample Recovery recommendations on page 11 for details of suggested dilutions. **If the end user chooses to use a different buffer as the Sample Diluent, it is up to the end user to determine if the buffer works in the assay, the appropriate dilution of samples, and assay validation.**

Samples in Tissue Culture Media can also be read in the assay provided the standards have been diluted into the Tissue Culture Media. There will be a small change in the standard curve associated with running the standards and samples in media. Media containing ferrous salts should be avoided as they will interfere with sensitive detections. The presence of pH indicator will not affect relative assay detection as long as standards and samples are in the same media.

PROCEDURAL NOTES

1. Do not mix components from different lot numbers or use reagents beyond the expiration date.
2. Allow all reagents, with the exception of the Hydrogen Peroxide Color Reagent, to warm to room temperature for at least 30 minutes before opening.
3. The Hydrogen Peroxide Color Reagent must be kept at 4°C during use.
4. Standards can be made up in either glass or plastic tubes.
5. Pre-rinse the pipet tip with the reagent, use fresh pipet tips for each sample, standard and reagent.
6. Pipet standards and samples to the bottom of the wells.
7. Add the reagent to the side of the well to avoid contamination.

REAGENT PREPARATION

Hydrogen Peroxide Standard

Allow the 100,000ng/mL Hydrogen Peroxide standard solution to warm to room temperature. Label six 12 x 75 mm tubes #1 through #6.

Pipet 966 μ l of sample diluent (buffer or Tissue Culture Media) into tube #1. Pipet 500 μ l of Diluent into tubes #2 - #6.

Add 34 μ l of the 100,000ng/mL standard to tube #1. Vortex thoroughly. Add 500 μ l of tube #1 to tube #2 and vortex thoroughly. Add 500 μ l of tube #2 to tube #3 and vortex. Continue this for tubes #4 through #6.

The concentration of Hydrogen Peroxide in tubes #1 through #6 will be 3,400, 1,700, 850, 425, 212.5 and 106.25ng/mL respectively. This converts to 100 μ M, 50 μ M, 25 μ M, 12.5 μ M, 6.25 μ M and 3.125 μ M respectively. See Hydrogen Peroxide Assay Layout Sheet for dilution details.

ASSAY PROCEDURE

Bring all reagents to room temperature for at least 30 minutes prior to opening. All standards and samples should be run in duplicate.

1. Determine the number of wells to be used. Cover unused wells tightly with a plate sealer. **DO NOT REUSE WELLS!**
2. Pipet 50 μ l of sample diluent (buffer or Tissue Culture Media) into duplicate Blank (Zero Standard) wells.
3. Pipet 50 μ l of Standards #1 through #6 into duplicate wells.
4. Pipet 50 μ l of Samples into duplicate wells.
5. Pipet 100 μ l of Color Reagent into the Blank, Standards and Sample wells.
6. Mix well by shaking or tapping the side of the plate for 10 seconds.
7. Incubate for 30 minutes at room temperature.
8. Blank the plate reader against the blank wells, read the optical density between 540 and 570nm, preferably at 550nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the blank wells from all readings.

CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of Hydrogen Peroxide in the samples. We recommend that the data be handled by an immunoassay software package utilizing a four parameter logistic curve fitting program (such as AssayBlaster™, catalog number ADI-28-0002). Such software is often supplied by plate reader manufacturers. If this type of data reduction software is not readily available, the concentration of Hydrogen Peroxide can be calculated as follows:

1. Calculate the average net Optical Density (OD) for each standard and sample by subtracting the average Blank OD from the average OD for each standard and sample:

$$\text{Average Net OD} = \text{Average OD} - \text{Average Blank OD}$$

2. Plot the Average Net OD for each Standard versus Hydrogen Peroxide Concentration. Approximate a straight line through the points. The concentration of Hydrogen Peroxide in the unknowns can be determined by interpolation.

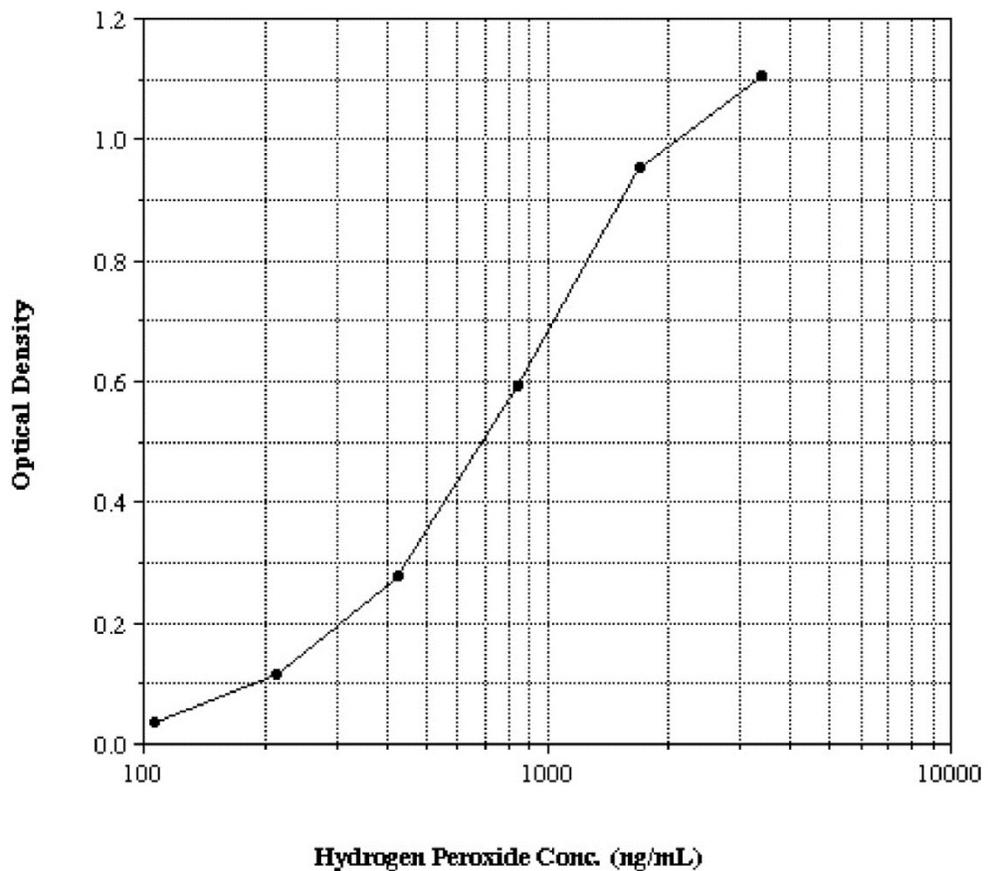
TYPICAL RESULTS

The results shown below are for illustration only and should not be used to calculate results.

Sample	Average OD	Net OD	Hydrogen Peroxide	
			(ng/mL)	μM
Blank	0.306			
S1	1.412	1.106	3,400	100
S2	1.262	0.956	1,700	50
S3	0.899	0.593	850	25
S4	0.585	0.279	425	12.5
S5	0.421	0.115	212.5	6.25
S6	0.341	0.035	106.25	3.125
Unknown1	0.838	0.532	767.6	22.58
Unknown2	1.200	0.894	1,555.1	45.74

TYPICAL STANDARD CURVES

Typical standard curves are shown below. These curves must not be used to calculate hydrogen peroxide concentrations; each user must run a standard curve for each assay.



PERFORMANCE CHARACTERISTICS

The following parameters for this kit were determined using the guidelines listed in the National Committee for Clinical Laboratory Standards (NCCLS) Evaluation Protocols¹⁴.

Sensitivity

Hydrogen Peroxide sensitivity in 50mM Phosphate, pH 6.0 was calculated by determining the average OD bound for sixteen (16) wells run as the zero standard, and comparing to the average OD for sixteen (16) wells run with Standard #6. The detection limit was determined as the concentration of Hydrogen Peroxide measured at two (2) standard deviations from the zero along the standard curve.

OD for Zero Standard =	0.277 ± 0.020 (7.3%)
OD for Standard #6 =	0.362 ± 0.051 (14.0%)
Delta OD (106.25-0ng/mL) = 0.362 - 0.277 =	0.085
2 SD's of Zero Standard =	0.041
Sensitivity = $\frac{0.041}{0.085} \times 106.25\text{ng/mL}$ =	51.25ng/mL

Linearity

A sample containing 2,880ng/mL Hydrogen Peroxide was serially diluted 5 times 1:2 in the recommended sample diluent and measured in the assay. The data was plotted graphically as actual Hydrogen Peroxide concentration versus measured Hydrogen Peroxide concentration. The line obtained had a slope of 0.9342 with a correlation coefficient of 0.9983.

Precision

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of Hydrogen Peroxide and running these samples multiple times (n=16) in the same assay. Inter-assay precision was determined by measuring three samples with low, medium and high concentrations of Hydrogen Peroxide in multiple assays (n=8). The precision numbers listed below represent the percent coefficient of variation for the concentrations of Hydrogen Peroxide determined in these assays as calculated by a curve fitting program.

	Hydrogen Peroxide (ng/mL)	Intra-assay %CV	Inter-assay %CV
Low	314.2		10.0
Medium	772.1		3.0
High	1,606		4.0
Low	326.5	2.2	
Medium	768.2	1.7	
High	1,732	5.7	

SAMPLE RECOVERIES

Please refer to pages **4-5** for Sample Handling recommendations and Standard preparation.

Hydrogen Peroxide concentrations were measured in horse heparinized plasma, human serum, human urine and tissue culture media. Hydrogen Peroxide was spiked into the undiluted samples of these matrices which were then diluted with the appropriate diluent and assayed in the kit. The following results were obtained:

Sample	% Recovery*	Recommended Dilution*
Horse heparin Plasma	94.5	>1:64
Human Serum	94.7	>1:64
Human Urine	90.6	>1:64
Tissue Culture Media	105.6	None

* See instructions in section titled "Sample Handling" for details.

REFERENCES

1. Davies KJ. IUBMB Life (1999) 48(1):41-7.
2. Zhang J et al. Antioxid. Redox Signal (2001) Jun;3(3):493-504.
3. Li, N & M. Karin Is NF- κ B the sensor of oxidative stress? FASEB J. 1999 13:1137-43.
4. Kim DK, et al. (2001) J. Cell Sci.. 114:4329-34.
5. Emelyanov A., et al. (Chest (2001) 120:1136-9.
6. Uesuge M., et al. J. Immunol. (2000) 165:6532-7.
7. Okuda M. et al. Artheroscler. Thromb. Vasc. Biol. (2001) Sep;21(9):1483-7.
8. Peiro C. et al. Br. J. Pharmacol. (2001) Aug;133(7):967-74.
9. Mody N. et al. Free Radic. Biol. Med. (2001) Aug;31(4):509-19.
10. Halliwell B. Drugs Aging (2001) 18(9):685-716.
11. Sanji E. et al. Biochem. Biophys. Res. Commun. (2001) Oct 5;287(4):1003-8.
12. Wentworth Jr. P. et al. Science (2001) 293:1806-11.
13. Wentworth, AD. Et al. PNAS (2000) 97(20):10930-5.
14. National Committee for Clinical Laboratory Standards Evaluation Protocols, SC1, 1989,NCCLS, Villanova, PA, 19085.



Product Manual

NOTES



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:

www.enzolifesciences.com