



Hydrogen peroxide chemiluminescent detection kit

Catalog No. ADI-907-012

96 Determination Kit

Table of Contents

Description	Page	2
Introduction		2
Precautions		2
Materials Supplied		3
Storage		3
Materials Needed but Not Supplied		3
Sample Handling		4
Procedural Notes		4
Reagent Preparation		5
Luminescent Assay Procedure		6
Calculation of Results		7
Typical Results		7
Typical Standard Curves		8
Performance Characteristics		9
References		11
Limited Warranty		12

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

Description

The Hydrogen peroxide chemiluminescent detection kit is a complete kit for the quantitative determination of Hydrogen Peroxide in colorless tissue culture media and buffers. Please read the complete kit insert before performing this assay.

Introduction

Hydrogen Peroxide (H_2O_2) 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 H_2O_2 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. The microtiter plate supplied in the kit has been selected for its low luminescent background and excellent reproducibility.
3. 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 Sample Handling, page 4.

Materials Supplied

1. **White Microtiter Plate, 1 each, Catalog No. 80-0953**
Break apart microtiter plate. Store at room temperature in the plastic ziploc bag provided to prevent contamination.
2. **Hydrogen Peroxide Standard, 0.5 mL, Catalog No. 80-0941**
A solution of Hydrogen Peroxide at 100,000 ng/mL in water with preservatives.
3. **Hydrogen Peroxide Substrate, 6 mL, Catalog No. 80-0939**
A solution of chemiluminescent substrate in aqueous buffers. **Substrate is sensitive to light. Store in the dark. Protect from ambient room light.**
4. **Hydrogen Peroxide Trigger Solution, 15 mL, Catalog No. 80-0940**
A solution of Bovine Hemoglobin containing preservatives.
5. **Hydrogen Peroxide Assay Layout Sheet, 1 each, Catalog No. 30-0165**

Storage

All components of this kit are stable at 4°C until the kit's expiration date.

Materials Needed but Not Supplied

1. Precision pipets for volumes between 25 μ L and 1,000 μ L.
2. Repeater pipet for dispensing 50 μ L.
3. Plate Luminometer with one injector capable of flash luminescence. For help with suitable instrumentation please contact the Technical Services department.
4. Graph paper for plotting the standard curve.

Sample Handling

The Hydrogen peroxide chemiluminescent detection kit is compatible with samples in Tissue Culture Media and buffer. Samples in these matrices can be read directly from the standard curve, provided the standards have been diluted in the same diluent as the samples. A tissue culture media **without** phenol red or other color indicators is recommended for use in this assay. **Phenol red has been shown to interfere.**

This assay is not designed for samples which could potentially contain Hemoglobin. The hydrogen peroxide in the standard or sample reacts with Hemoglobin, which in this system is used as the trigger. The product of this reaction then reacts with the provided substrate, yielding light as a final product. The problem with biological samples containing Hemoglobin is it will trigger the Substrate before the plate is placed in the plate reader and the reaction is measured.

An alternative method for biological samples would be to directly pipet the standards or samples and Hemoglobin into the wells and use the Substrate as the trigger in the luminometer. However, after using this method, the trigger lines used will have to be designated for this Substrate only and will have to be replaced for other triggers. This alternative method is offered only as a suggestion.

Procedural Notes

1. Do not mix components from different lot numbers or use reagents beyond the expiration date.
2. Allow all reagents to warm to room temperature for at least 30 minutes before opening.
3. Standards can be made up in either glass or plastic tubes.
4. Pre-rinse the pipet tip with the reagent, use fresh pipet tips for each sample, standard and reagent.
5. Pipet standards and samples to the bottom of the wells.
6. Add the reagents to the side of the well to avoid contamination.
7. This kit uses break-apart microtiter strips, which allow the user to measure as many samples as desired. Unused wells can be kept at room temperature in the plastic ziploc bag provided. The wells should be used in the frame provided.
8. **This assay uses a Luminescent measurement of Hydrogen Peroxide concentration. The luminescent signal is typically represented as Relative Light Units (RLU). Different luminometers will display different RLU readings. Please see the luminometer instruction manual for details.**

Reagent Preparation

1. Hydrogen Peroxide Standard(s)

In buffer:

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

Pipet 975 μ L of buffer into tube #1. Pipet 750 μ L of buffer into tubes #2 - #6.

Add 25 μ L of the 100,000 ng/mL standard to tube #1. Vortex thoroughly. Add 250 μ L of tube #1 to tube #2 and vortex thoroughly. Add 250 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 2,500, 625, 156.25, 39.06, 9.77 and 2.44 ng/mL respectively. See Hydrogen Peroxide Assay Layout Sheet for dilution details.

In Tissue Culture Media:

Allow the 100,000 ng/mL Hydrogen Peroxide Standard Solution to warm to room temperature. Label five 12 x 75 mm glass tubes #1 through #5.

Pipet 900 μ L of tissue culture media into tube #1. Pipet 750 μ L of tissue culture media into tubes 2 - #5.

Add 100 μ L of the 100,000 ng/mL standard to tube #1. Vortex thoroughly. Add 250 μ L of tube #1 to tube #2 and vortex thoroughly. Continue this for tubes #3 through #5.

The concentration of Hydrogen Peroxide in tubes #1 through #5 will be 10,000, 2,500, 625, 156.25 and 39.06 ng/mL respectively. See Hydrogen Peroxide Assay Layout Sheet for dilution details.

Assay Procedure

All samples should be run in duplicate.

All samples should be allowed to warm to room temperature for at least 30 minutes prior to use.

1. Determine the number of wells to be used and put any remaining wells back into plastic ziploc bag. Store unused wells at room temperature.
2. Pipet 50 μ L of standard diluent (buffer or tissue culture media) into duplicate Blank (Zero Standard) wells.
3. Pipet 50 μ L of Standards #1 through #6 for buffer, or #1 through #5 for tissue culture media, into duplicate wells.
4. Pipet 50 μ L of Samples into duplicate wells.
5. Pipet 50 μ L of Substrate into the Blank, Standards and Sample wells.
6. Mix well by shaking or tapping the side of the plate for 10 seconds.
7. Place microtiter plate in luminometer for chemiluminescent measurement.
8. Inject 50 μ L of Hydrogen Peroxide Trigger into each well. Immediately read in luminometer for 5 seconds.
9. Determine integrated light output for the 5 second read time in Relative Light Units (RLU).

For Step 8: Program your luminometer to inject with zero delay between injection and light detection. Read time should be set at 5 seconds.

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 a software package utilizing a 4 parameter logistic curve fitting program. 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 Relative Light Units (RLU) for each standard and sample by subtracting the average Blank RLU from the average RLU for each standard and sample:

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

2. Plot the Average Net RLU 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

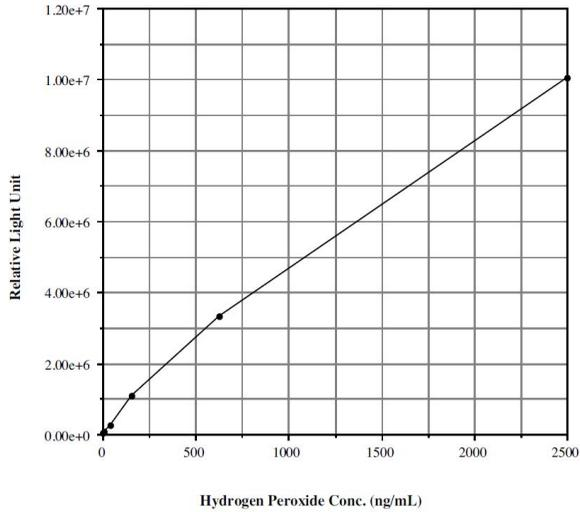
The results shown below are for illustration only and **should not** be used to calculate results from another assay.

	<u>BUFFER</u>			<u>TISSUE CULTURE MEDIA</u>		
<u>Sample</u>	<u>Average RLU</u>	<u>Net RLU</u>	<u>Hydrogen Peroxide ng/mL</u>	<u>Average RLU</u>	<u>Net RLU</u>	<u>Hydrogen Peroxide ng/mL</u>
Blank	(129,892)			(146,118)		
S1	10,160,143	10,030,251	2,500	2,291,888	2,145,770	10,000
S2	3,462,480	3,332,588	625	598,190	452,072	2,500
S3	1,228,303	1,098,411	156.25	247,691	101,573	625
S4	396,837	266,945	39.06	165,902	19,784	156.25
S5	201,917	72,025	9.77	150,207	4,089	39.06
S6	157,039	27,147	2.44			

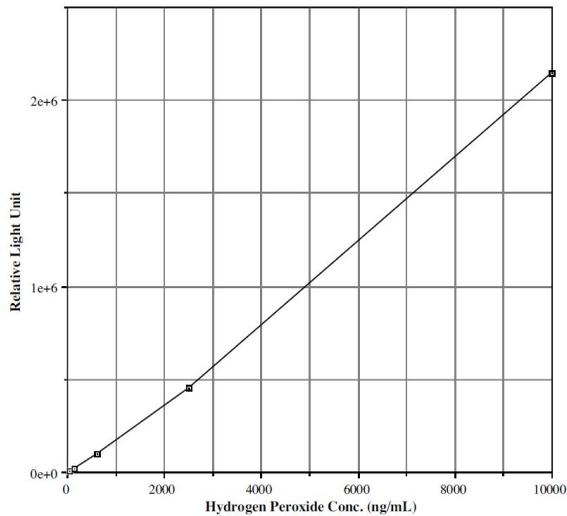
Typical Standard Curves

Typical standard curves are shown below. The curves **must not** be used to calculate Hydrogen Peroxide concentrations; each user must run a standard curve for each assay.

Standard Curve in Buffer



Standard Curve in Tissue Culture Media



Performance Characteristics

The following parameters for this kit were determined using the guidelines listed in the National Committee for Clinical Laboratory Standards (NCCLS) Evaluation Protocols.¹⁴

Sensitivity

Hydrogen Peroxide sensitivity was calculated by determining the average RLU bound for sixteen (16) wells run as the zero standard, and comparing to the average RLU for sixteen (16) wells run with Standard #6 for Buffer or Standard #5 for Tissue Culture Media. The detection limit was determined as the concentration of Hydrogen Peroxide measured at two (2) standard deviations from the zero along the standard curve.

Buffer Sensitivity

RLU for Zero Standard = 129,168 ± 9,374 (7.3%)

RLU for Standard #6 = 148,271 ± 5,254 (3.5%)

Delta RLU (2.44-0 ng/mL) = 148,271-129,168 = 19,102

2 SD's of S0 = 18,748

Sensitivity = $\frac{18,748}{19,102} \times 2.44 \text{ ng/mL} = \mathbf{2.39 \text{ ng/mL}}$

Tissue Culture Media

RLU for Zero Standard = 143,039 ± 2,949 (2.1%)

RLU for Standard #6 = 162,409 ± 3,513 (2.2%)

Delta RLU (39.1-0 ng/mL) = 162,409-143,039 = 19,370

2 SD's of S0 = 5,898

Sensitivity = $\frac{5,898}{19,370} \times 39.1 \text{ ng/mL} = \mathbf{11.91 \text{ ng/mL}}$

Linearity

Buffer

A sample containing 1,146 ng/mL Hydrogen Peroxide was serially diluted 7 times 1:2 in buffer 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 0.9841 with a correlation coefficient of 0.9998.

Tissue Culture Media

A sample containing 8,184 ng/mL Hydrogen Peroxide was serially diluted 6 times 1:2 in Tissue Culture Media 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 0.9441 with a correlation coefficient of 0.9997.

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.

	<u>BUFFER</u>			<u>TISSUE CULTURE MEDIA</u>		
	Hydrogen Peroxide (ng/mL)	Inter-assay (%CV)	Intra-assay (%CV)	Hydrogen Peroxide (ng/mL)	Inter-assay (%CV)	Intra-assay (%CV)
Low	27.33	7.8		33.1	8.9	
Medium	55.25	6.1				
High	1,135.85	8.4		880.4	3.9	
Low	26.49		9.6	483.04		3.4
Medium	51.57		6.7			
High	1,167.56		5.8	911.85		4.0

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? *FASEBJ.* 1999 13:1137-43. Review.
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) Aug165;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.



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 purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. 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

www.enzolifesciences.com
Enabling Discovery in Life Science®

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 NONINFRINGEMENT 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.

Global Headquarters

Enzo Life Sciences Inc.

10 Executive Blvd
Farmingdale, NY 11735
(p) 1-800-942-0430
(f) 1-631-694-7501
(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach
CH-4415 Lause / Switzerland
(p) +41/0 61 926 89 89
(f) +41/0 61 926 89 79
(e) info-ch@enzolifesciences.com

Please visit our website at www.enzolifesciences.com for additional contact information.

