

# Hydrogen Peroxide Fluorescent Detection Kit



Kit #9131, 2 plates  
**RESEARCH USE ONLY**

## 1. Introduction

Hydrogen peroxide was first described in 1818 by Louis Jacques Thénard. Today, hydrogen peroxide is industrially manufactured almost exclusively by the autoxidation of a 2-alkyl-9,10-dihydroxyanthracene to the corresponding 2-alkyl anthraquinone in the Riedl-Pfleiderer or anthraquinone process.

In biological systems, incomplete reduction of  $O_2$  during respiration produces superoxide anion ( $O_2^{\cdot-}$ ), which is spontaneously or enzymatically dismutated by superoxide dismutase to  $H_2O_2$ . Many cells produce low levels of  $O_2^{\cdot-}$  and  $H_2O_2$  in response to a variety of extracellular stimuli, such as cytokines (TGF- $\beta$ 1, TNF- $\alpha$ , and various interleukins), peptide growth factors (PDGF; EGF, VEGF, bFGF, and insulin), the agonists of heterotrimeric G protein-coupled receptors (GPCR) such as angiotensin II, thrombin, lysophosphatidic acid, sphingosine 1-phosphate, histamine, and bradykinin, and by shear stress<sup>1</sup>. The addition of exogenous  $H_2O_2$  or the intracellular production in response to receptor stimulation affects the function of various proteins, including protein kinases, protein phosphatases, transcription factors, phospholipases, ion channels, and G proteins. In 1894, Fenton<sup>2</sup> described the oxidation of tartaric acid by  $Fe^{2+}$  and  $H_2O_2$ .  $H_2O_2$  and  $O_2$  may participate in the production of singlet oxygen and peroxynitrite and the generation of these species may be concurrent with reactions involving iron, and under some circumstances, they might be important contributors to  $H_2O_2$  toxicity<sup>3,4</sup>.

A substantial portion of  $H_2O_2$  lethality involves DNA damage by oxidants generated from iron-mediated Fenton reactions<sup>5,6</sup>. Damage by Fenton oxidants occurs at the DNA bases or at the sugar residues. Sugar damage is initiated by hydrogen abstraction from one of the deoxyribose carbons, and the predominant consequence is eventual strand breakage and base release<sup>7,8</sup>.

ICT's Hydrogen Peroxide Fluorescent Detection Kit is designed to quantitatively measure  $H_2O_2$  in a variety of samples. This kit is validated for use in fresh urine, buffers, and tissue culture media (TCM) samples. It is species independent. Please read the complete kit insert before performing this assay. A hydrogen peroxide standard is provided to generate a standard curve for the assay. All samples should be read off the standard curve. Samples are mixed with the Fluorescent  $H_2O_2$  Detection Substrate and the reaction is initiated by addition of horseradish peroxidase. The reaction is incubated at room temperature for 15 minutes. HRP is oxidized by hydrogen peroxide present in the sample. Oxidized HRP then reacts with the substrate to convert the colorless substrate into the fluorescent form. The fluorescent product is read at 590 nm with excitation at 570 nm. Increasing levels of  $H_2O_2$  cause a linear increase in fluorescent product. This kit is for research use only and is not for use in diagnostic procedures.

*Learn more about all of ICT's products at  
[www.immunochemistry.com](http://www.immunochemistry.com) or call 1-800-829-3194.*

## 2. Kit Contents

- 2 black half-area polystyrene 96-well microtiter plates #268
- 1 vial of Hydrogen Peroxide Standard (220  $\mu$ L) #6603: Hydrogen Peroxide at 100  $\mu$ M in a special stabilizing solution (Section 12).
- 1 bottle of Assay Buffer Concentrate (25 mL) #6605: A 5X buffer concentrate containing detergents and stabilizers (Section 10).
- 1 vial of Fluorescent  $H_2O_2$  Detection Substrate (5 mL) #6606: A solution of the substrate in a special stabilizing buffer.
- 1 vial of Horseradish Peroxidase Concentrate (60  $\mu$ L) #6607: A 100X concentrated solution of HRP in a special stabilizing solution (Section 11).

## 3. Required Materials

- Repeater pipet with disposable tips capable of accurately dispensing 25  $\mu$ L.

## 4. Storage

All components of this kit should be stored at 4°C until the expiration date of the kit.

## 5. MSDS

Material safety data sheets are available online at [www.immunochemistry.com](http://www.immunochemistry.com) or by calling 1-800-829-3194 or 952-888-8788.

## 6. Precautions

As with all such products, this kit should only be used by qualified personnel who have had laboratory safety instruction. The complete insert should be read and understood before attempting to use the product. The supplied hydrogen peroxide standard consists of a very dilute  $H_2O_2$  solution.

## 7. Detection Equipment

The assay can be analyzed with a fluorescence plate reader:

- 96-well plate reader capable of reading fluorescence at 580-590 nm with excitation at 570-580 nm. Set plate parameters for a 96-well Corning® Costar 3694 plate.
- Software for converting fluorescent intensity readings from the plate reader and carrying out four parameter logistic curve (4PLC) fitting. Contact your plate reader manufacturer for details.

## 8. Sample Types and Preparation

Samples that need to be stored after collection should be stored at -70°C or lower, preferably after being frozen in liquid nitrogen. Urine samples can be used after being diluted  $\geq$  1:10. This assay has been validated for buffer and media samples.

	<b>1/2 Plate</b>	<b>1 Plate</b>	<b>2 Plates</b>
<b>Horseradish Peroxidase Concentrate</b>	15 $\mu$ L	30 $\mu$ L	55 $\mu$ L
<b>Assay Buffer</b>	1.485 mL	2.97 mL	5.445 mL
<b>Total Volume</b>	1.5 mL	3 mL	5.5 mL

	<b>Std 1</b>	<b>Std 2</b>	<b>Std 3</b>	<b>Std 4</b>	<b>Std 5</b>	<b>Std 6</b>	<b>Std 7</b>
<b>Assay Buffer (<math>\mu</math>L)</b>	450	200	200	200	200	200	200
<b>Addition</b>	Stock	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6
<b>Vol of Addition (<math>\mu</math>L)</b>	50	200	200	200	200	200	200
<b>Final Conc (<math>\mu</math>M)</b>	10	5	2.5	1.25	0.625	0.313	0.1569

### 9. Sample Preparation

Dilute samples  $\geq$  1:10 with Assay Buffer prior to running in the assay.

### 10. Assay Buffer Preparation

Dilute Assay Buffer Concentrate (#6605) 1:5 by adding one part of the concentrate to four parts of deionized water. Diluted Assay Buffer is stable at 4°C for 3 months.

### 11. Horseradish Peroxidase (HRP) Preparation

Dilute Horseradish Peroxidase Concentrate (#6607) 1:100 with Assay Buffer. See Table 1: HRP Dilution; for example if using 1 plate:

1. Measure 2.97 mL Assay Buffer.
2. Add 30  $\mu$ L Horseradish Peroxidase Concentrate.
3. Mix. The total volume is 3 mL for 1 plate.

### 12. Standard Preparation

1. Hydrogen peroxide standards are prepared by labeling seven tubes as #1 through #7.
2. Briefly vortex to mix the vial of H<sub>2</sub>O<sub>2</sub> Standard (#6603).
3. Pipet 450  $\mu$ L of Assay Buffer into tube #1 and 200  $\mu$ L into tubes #2 to #7.
4. Carefully add 50  $\mu$ L of the H<sub>2</sub>O<sub>2</sub> Standard to tube #1 and vortex completely.
5. Take 200  $\mu$ L of the solution in tube #1 and add it to tube #2 and vortex completely.
6. Repeat this for tubes #3 through #7.
7. The concentration of H<sub>2</sub>O<sub>2</sub> in tubes 1 through 7 will be 10, 5, 2.5, 1.25, 0.625, 0.313 and 0.156  $\mu$ M (See Table 2: Standard Preparation).

- Use all Standards within 2 hours of preparation.

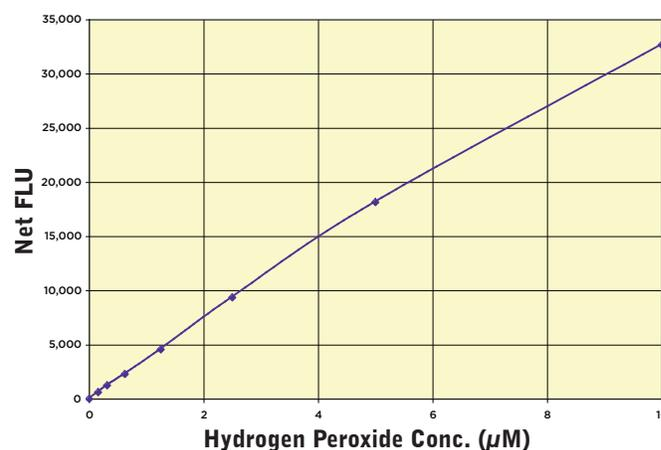
**Table 3: Typical Data**

Always run your own standard curves for calculation of results. Do not use this data. Conversion Factor: 100 nM of hydrogen peroxide is equivalent to 3.4 ng/mL.

<b>Sample</b>	<b>Mean FLU</b>	<b>Net FLU</b>	<b>H<sub>2</sub>O<sub>2</sub> Conc. (<math>\mu</math>M)</b>
<b>Zero</b>	3,782	0	0
<b>Standard 1</b>	36,417	32,635	10
<b>Standard 2</b>	21,919	18,137	5
<b>Standard 3</b>	13,134	9,352	2.5
<b>Standard 4</b>	8,333	4,551	1.25
<b>Standard 5</b>	6,072	2,290	0.625
<b>Standard 6</b>	5,031	1,249	0.313
<b>Standard 7</b>	4,398	616	0.156
<b>Sample 1</b>	6,578	2,796	0.76
<b>Sample 2</b>	24,680	20,898	5.85

**Figure 1: Typical Standard Curve**

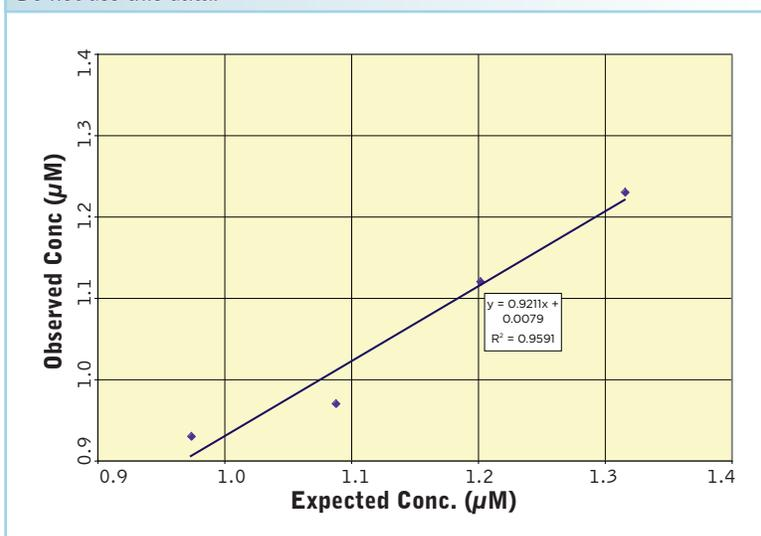
Always run your own standard curves for calculation of results. Do not use this data.



High H <sub>2</sub> O <sub>2</sub> RPMI	Low H <sub>2</sub> O <sub>2</sub> RPMI	Observed Conc. (μM)	Expected Conc. (μM)	% Recovery
80%	20%	1.23	1.32	93.5
60%	40%	1.12	1.20	93.2
40%	60%	0.97	1.09	89.2
20%	80%	0.93	0.97	95.5
<b>Mean Recovery</b>				<b>92.8%</b>

**Figure 2: Sample Linearity Curve**

Always run your own standard curves for calculation of results.  
Do not use this data.



### 13. Assay Protocol

1. Use the plate layout sheet on the back page to aid in proper sample and standard identification. Set plate parameters for a 96-well Corning Costar 3694 plate.
2. Pipet 50 μL of samples or appropriate standards into duplicate wells in the plate.
3. Pipet 50 μL of Assay Buffer into duplicate wells as the zero standard.
4. Add 25 μL of Fluorescent H<sub>2</sub>O<sub>2</sub> Detection Substrate (#6606) to each well using a repeater pipet.
5. Initiate the reaction by adding 25 μL of the HRP Preparation (Section 11) to each well using a repeater pipet. Replace tips each time to minimize sample carry over.
6. Incubate at room temperature for 15 minutes.
7. Read the fluorescent emission at 585 ± 5 nm with excitation at 575 ± 5 nm. Please contact your plate reader manufacturer for suitable filter sets.

### 14. Calculation of Results

1. Set up fluorescence plate reader software to subtract mean of the zero 0-well FLU values from all standard and sample fluorescence readings. This 0-well FLU subtraction step can be performed after the plate reader has completed the fluorescence scan of the plate.
2. Manually or via plate reader software, calculate the average FLU reading for each of the duplicate standards and samples.
3. Create a standard curve using software-derived linear regression analysis. Select the four-parameter logistic curve (4PLC) fitting option for this step (see Figure 1: Typical Standard Curve).

4. Multiply curve derived sample concentration values by initial sample dilution factor to obtain the H<sub>2</sub>O<sub>2</sub> concentration present in neat samples.

### 15. Validation Data: Sensitivity

Sensitivity was calculated by comparing the FLUs for twenty wells run for each of the zero and standard #7 (low standard). The theoretical detection limit in Assay Buffer was determined at two (2) standard deviations from the zero along the standard curve.

Sensitivity was determined as 0.038 μM. This is equivalent to 1.9 pmol (64.6 pg) H<sub>2</sub>O<sub>2</sub> per well.

### 16. Validation Data: Limit of Detection

The limit of detection in a human sample was determined in a similar manner by comparing the FLUs for twenty wells run for each of the zero and a low concentration human sample.

The Limit of Detection was determined as 0.052 μM. This is equivalent to 2.6 pmol (88.4 pg) H<sub>2</sub>O<sub>2</sub> per well.

### 17. Linearity

Linearity was determined by taking two RPMI-1640 media samples with known H<sub>2</sub>O<sub>2</sub> concentrations and mixing them in the ratios given; see Table 4: Sample Linearity Data. The measured concentrations were compared to the expected values based on the ratios used. Figure 2 illustrates a linear plot of observed versus expected concentration values.

**Table 5: Intra Assay Precision**

Three buffer samples were run in replicates of 20 in an assay. The mean and precision of the calculated concentrations were:

Sample	H <sub>2</sub> O <sub>2</sub> Conc. (μM)	%CV
1	6.27	3.6
2	3.21	3.8
3	0.98	5.7

**Table 6: Inter Assay Precision**

Three buffer samples were run in duplicates in fourteen assays run over multiple days by three operators. The mean and precision of the calculated concentrations were:

Sample	H <sub>2</sub> O <sub>2</sub> Conc. (μM)	%CV
1	5.86	4.3
2	3.00	7.0
3	0.88	12.1

**18. References**

1. Rhee, S.G., et al. Hydrogen peroxide: A key messenger that modulates protein phosphorylation through cysteine oxidation. *Science's stke*. (2000). Available at: <http://stke.sciencemag.org/cgi/content/abstract/sigtrans;2000/53/pe1>
2. Fenton, H.J.H. *J. Chem. Soc. (Lond.)* 65, 899–910. (1894).
3. Sies, H. *Mutat. Res.* 299, 183–191. (1993).
4. Squadrito, G.L. and Pryor, W.A. The formation of peroxynitrite *in vivo* from nitric oxide and superoxide. *Chem. Biol. Interact.* 96, 203–206. (1995).
5. Imlay, J.A. and Linn, S. DNA damage and oxygen radical toxicity. *Science* 240, 1302–1309. (1988).
6. Mello-Filho, A.C. and Meneghini, R. Iron is the intracellular metal involved in the production of DNA damage by oxygen radicals. *Mutat. Res.*, 251, 109–113. (1991).
7. von Sonntag, C. In: *The Chemical Basis of Radiation Biology*. pp. 238–249, Taylor and Francis, New York. (1987).
8. Henle, E.S. et al. DNA strand breakage is correlated with unaltered base release after gamma irradiation. *Radiat. Res.* 143, 144–150. (1995).

	1	2	3	4	5	6	7	8	9	10	11	12
A												
B												
C												
D												
E												
F												
G												
H												

*Thank you for using this kit. If you have any questions or would like to share your data, please contact us at 1-800-829-3194 or 952-888-8788, or send an email to [help@immunochemistry.com](mailto:help@immunochemistry.com).*

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



ImmunoChemistry Technologies, LLC  
9401 James Avenue South #155  
Bloomington, MN 55431 USA

toll-free: 1-800-829-3194  
local: 952-888-8788  
fax: 952-888-8988  
[www.immunochemistry.com](http://www.immunochemistry.com)