

## Hydrogen Peroxide Colorimetric/Fluorometric Assay Kit

(Catalog #K265-200; 200 reactions; Store kit at -20°C)

### I. Introduction:

Hydrogen Peroxide is a reactive oxygen metabolic byproduct that serves as a key regulator for a number of oxidative stress-related states. Functioning through NF-κB and other factors, hydroperoxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis, diabetic vasculopathy, osteoporosis, neuro-degenerative diseases, Down's syndrome and immune system diseases. **BioVision's Hydrogen Peroxide Assay Kit** is a highly sensitive, simple, direct and HTS-ready colorimetric and fluorometric assay for measuring H<sub>2</sub>O<sub>2</sub> in biological samples. In the presence of Horse Radish Peroxidase (HRP), the OxiRed Probe reacts with H<sub>2</sub>O<sub>2</sub> to produce product with color ( $\lambda_{max}$  = 570 nm) and red-fluorescent (Ex/Em = 535/587 nm). The kit can perform 200 reactions by fluorometric method or 100 reactions by colorimetric method. The detection limit can be as low as 2 pmol/well (or 40 μM) of H<sub>2</sub>O<sub>2</sub> in the sensitive fluorometric assay.

### II. Kit Contents:

Components	K265-200	Cap Code	Part Number
H <sub>2</sub> O <sub>2</sub> Assay Buffer	25 ml	WM	K265-200-1
OxiRed™ Probe (in anhydrous DMSO)	0.2 ml	Red	K265-200-2A
HRP	1 vial	Green	K265-200-4
H <sub>2</sub> O <sub>2</sub> Standard (0.88 M)	0.1 ml	Yellow	K265-200-5

### III. Storage and Handling:

Warm the assay buffer to room temperature (RT) before use. Briefly centrifuge vials prior to opening. Read the entire protocol before performing the assay.

### IV. Reagent Reconstitution and General Consideration:

**OxiRed™ Probe:** Ready to use as supplied. Briefly warm at 37°C to melt the frozen DMSO.

The OxiRed™ Probe solution is stable for 1 week at 4°C and 1 month at -20°C.

**HRP:** Dissolve in 220 μl H<sub>2</sub>O<sub>2</sub> Assay Buffer, pipetting up and down. The HRP solution is stable for 1 week at 4°C and 1 month at -20°C.

### V. Hydrogen Peroxide Assay:

#### 1. Sample Preparation:

Collect the cell culture supernatant, serum, plasma, urine and other biological fluids (contains 0.8 - 6 μM H<sub>2</sub>O<sub>2</sub>). Centrifuge for 15 min at 1000 x g within 30 min of collection. Remove the particulate pellet. Samples, especially those such as culture medium, tissue lysate or plasma should be filtered through 10 kDa spin columns (BioVision, Cat# 1997-25) to remove all the proteins then kept at -80°C for storage. It is recommended to assay the sample(s) immediately or aliquot and store at -80°C. Avoid repeated freeze-thaw cycles. Add 2 - 50 μl samples into each well. Bring the volume to 50 μl with H<sub>2</sub>O<sub>2</sub> Assay Buffer.

#### 2. H<sub>2</sub>O<sub>2</sub> Standard Curve:

**For the Colorimetric Assay:** Dilute 10 μl 0.88 M H<sub>2</sub>O<sub>2</sub> Standard into 870 μl dH<sub>2</sub>O to generate a 10 mM H<sub>2</sub>O<sub>2</sub> Standard. Then dilute 10 μl of the 10 mM H<sub>2</sub>O<sub>2</sub> Standard into 990 μl dH<sub>2</sub>O to generate a 0.1 mM H<sub>2</sub>O<sub>2</sub> Standard. Add 0, 10, 20, 30, 40, 50 μl of the 0.1 mM H<sub>2</sub>O<sub>2</sub> Standard into 96-well plate in duplicate to generate 0, 1, 2, 3, 4, 5 nmol/well H<sub>2</sub>O<sub>2</sub> Standard. Adjust the volume to 50 μl/well using H<sub>2</sub>O.

**For the Fluorometric Assay:** Dilute 100 μl of the 0.1 mM H<sub>2</sub>O<sub>2</sub> Standard into 900 μl dH<sub>2</sub>O to generate a 10 μM H<sub>2</sub>O<sub>2</sub> Standard. Add 0, 10, 20, 30, 40, 50 μl of the 10 μM H<sub>2</sub>O<sub>2</sub> Standard into a 96-well plate in duplicate to generate 0, 0.1, 0.2, 0.3, 0.4, 0.5 nmol/well H<sub>2</sub>O<sub>2</sub> Standard. Adjust the volume to 50 μl/well using H<sub>2</sub>O.

**3. Reaction Mix:** Mix enough reagents for the number of assays to be performed. For each well, prepare a total 50 μl Reaction Mix:

#### Colorimetric Assay

46 μl Assay Buffer  
2 μl OxiRed™ Probe solution  
2 μl HRP solution

#### Fluorometric Assay

48 μl Assay Buffer  
1 μl OxiRed™ Probe solution  
1 μl HRP solution

Add 50 μl of the Reaction Mix to each test samples and H<sub>2</sub>O<sub>2</sub> Standard wells. Mix well. Incubate at RT for 10 min.

\*For a more sensitive assay, you can dilute the Standard 10 fold further, decrease the OxiRed™ amount to 0.2 μl and HRP amount to 0.4 μl per well,. This will decrease the fluorescence background and detects as low as 2 pmol/well (or 40 μM) H<sub>2</sub>O<sub>2</sub>.

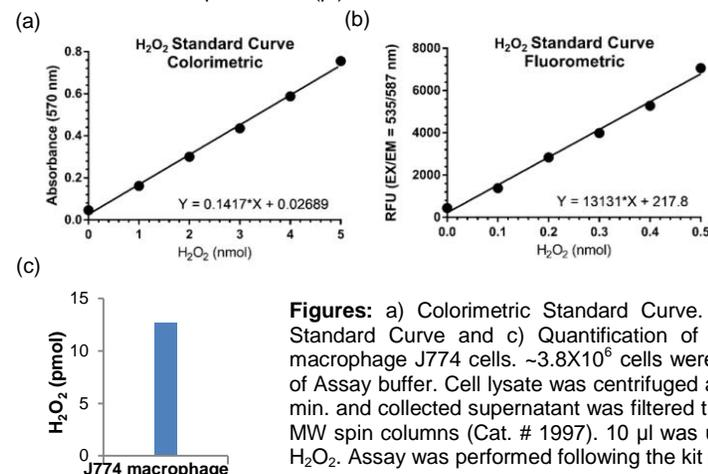
**4. Measure OD (570 nm) or fluorescence (Ex/Em = 535/587 nm) in a micro-plate reader.**

**5. Calculation:** Correct background by subtracting the value derived from the 0 nmol H<sub>2</sub>O<sub>2</sub> control from all sample and Standard readings (Note: The background reading can be significant and must be subtracted from sample readings). Plot the H<sub>2</sub>O<sub>2</sub> Standard Curve. Apply your sample readings to the Standard Curve. H<sub>2</sub>O<sub>2</sub> concentrations of the test samples can then be calculated,

$$C = Sa/Sv \text{ (pmol/}\mu\text{l or } \mu\text{M)},$$

Where: **Sa** is the sample amount from your Standard Curve (in pmol),

**Sv** is sample volume (μl).



**Figures:** a) Colorimetric Standard Curve. b) Fluorometric Standard Curve and c) Quantification of H<sub>2</sub>O<sub>2</sub> in murine macrophage J774 cells. ~3.8X10<sup>6</sup> cells were lysed in 380 μl of Assay buffer. Cell lysate was centrifuged at 10,000xg for 2 min. and collected supernatant was filtered through a 10 kDa MW spin columns (Cat. # 1997). 10 μl was used to measure H<sub>2</sub>O<sub>2</sub>. Assay was performed following the kit protocol.

### VI. RELATED PRODUCTS:

Glutathione Reductase Assay Kit  
Colorimetric Glutathione Detection Kit (GSH, GSSG and Total)  
GST Colorimetric Assay Kit  
Acid Phosphatase Assay Kit  
Phosphate Fluorescence Assay Kit  
NAD/NADH Quantification Kit

Glutathione Peroxidase Assay Kit  
ApoGSH Glutathione Detection Kit  
GST Fluorometric Assay Kit  
Triglyceride Assay Kit  
ADP/ATP Ratio Assay Kit  
Phosphate Colorimetric Assay Kit  
NADP/NADPH Quantitation Kit

**FOR RESEARCH USE ONLY! Not to be used on humans.**

## GENERAL TROUBLESHOOTING GUIDE:

Problems	Cause	Solution
Assay not working	<ul style="list-style-type: none"> <li>• Use of ice-cold assay buffer</li> <li>• Omission of a step in the protocol</li> <li>• Plate read at incorrect wavelength</li> <li>• Use of a different 96-well plate</li> </ul>	<ul style="list-style-type: none"> <li>• Assay buffer must be at room temperature</li> <li>• Refer and follow the data sheet precisely</li> <li>• Check the wavelength in the data sheet and the filter settings of the instrument</li> <li>• Fluorescence: Black plates ; Luminescence: White plates; Colorimeters: Clear plates</li> </ul>
Samples with erratic readings	<ul style="list-style-type: none"> <li>• Use of an incompatible sample type</li> <li>• Samples prepared in a different buffer</li> <li>• Samples were not deproteinized (if indicated in datasheet)</li> <li>• Cell/ tissue samples were not completely homogenized</li> <li>• Samples used after multiple free-thaw cycles</li> <li>• Presence of interfering substance in the sample</li> <li>• Use of old or inappropriately stored samples</li> </ul>	<ul style="list-style-type: none"> <li>• Refer data sheet for details about incompatible samples</li> <li>• Use the assay buffer provided in the kit or refer data sheet for instructions</li> <li>• Use the 10 kDa spin cut-off filter or PCA precipitation as indicated</li> <li>• Use Dounce homogenizer (increase the number of strokes); observe for lysis under microscope</li> <li>• Aliquot and freeze samples if needed to use multiple times</li> <li>• Troubleshoot if needed, deproteinize samples</li> <li>• Use fresh samples or store at correct temperatures till use</li> </ul>
Lower/ Higher readings in Samples and Standards	<ul style="list-style-type: none"> <li>• Improperly thawed components</li> <li>• Use of expired kit or improperly stored reagents</li> <li>• Allowing the reagents to sit for extended times on ice</li> <li>• Incorrect incubation times or temperatures</li> <li>• Incorrect volumes used</li> </ul>	<ul style="list-style-type: none"> <li>• Thaw all components completely and mix gently before use</li> <li>• Always check the expiry date and store the components appropriately</li> <li>• Always thaw and prepare fresh reaction mix before use</li> <li>• Refer datasheet &amp; verify correct incubation times and temperatures</li> <li>• Use calibrated pipettes and aliquot correctly</li> </ul>
Readings do not follow a linear pattern for Standard curve	<ul style="list-style-type: none"> <li>• Use of partially thawed components</li> <li>• Pipetting errors in the standard</li> <li>• Pipetting errors in the reaction mix</li> <li>• Air bubbles formed in well</li> <li>• Standard stock is at an incorrect concentration</li> <li>• Calculation errors</li> <li>• Substituting reagents from older kits/ lots</li> </ul>	<ul style="list-style-type: none"> <li>• Thaw and resuspend all components before preparing the reaction mix</li> <li>• Avoid pipetting small volumes</li> <li>• Prepare a master reaction mix whenever possible</li> <li>• Pipette gently against the wall of the tubes</li> <li>• Always refer the dilutions in the data sheet</li> <li>• Recheck calculations after referring the data sheet</li> <li>• Use fresh components from the same kit</li> </ul>
Unanticipated results	<ul style="list-style-type: none"> <li>• Measured at incorrect wavelength</li> <li>• Samples contain interfering substances</li> <li>• Use of incompatible sample type</li> <li>• Sample readings above/below the linear range</li> </ul>	<ul style="list-style-type: none"> <li>• Check the equipment and the filter setting</li> <li>• Troubleshoot if it interferes with the kit</li> <li>• Refer data sheet to check if sample is compatible with the kit or optimization is needed</li> <li>• Concentrate/ Dilute sample so as to be in the linear range</li> </ul>
<p><b>Note:</b> The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.</p>		