

# Catalase Activity Assay Kit

PromoKine

HTS-ready colorimetric & fluorometric assay for measuring catalase activity in biological samples

## Instruction Manual

| Catalog Number                                   | PK-CA577-K773   |            |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
|--|---|------------|----------|------------|-----------------------|-------|----|------------------------|--------|-----|-------------------|--------|-------|--|-------|--------|---------------|------|-------|---------------------------|------|------|
| Description                                      | <p>Catalase (EC 1.11.1.6) is a ubiquitous antioxidant enzyme that is present in nearly all living organisms. It functions to catalyze the decomposition of hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) to water and oxygen.</p> <p>PromoKine's Catalase Assay Kit provides a highly sensitive, simple, direct and HTS-ready assay for measuring Catalase activity in biological samples. In the assay, catalase first reacts with H<sub>2</sub>O<sub>2</sub> to produce water and oxygen, the unconverted H<sub>2</sub>O<sub>2</sub> reacts with OxiRed probe to produce a product, which can be measured at 570 nm (Colorimetric method) or at Ex/Em = 535/587 nm (fluorometric method). Catalase activity is reversely proportional to the signal. The kit can detect 1 μU or less of catalase activity in samples.</p>  |            |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Quantity   | 100 assays  |            |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Kit Components                                   | <table border="1"><thead><tr><th>Components</th><th>Quantity</th><th>Color Code</th></tr></thead><tbody><tr><td>Catalase Assay Buffer</td><td>25 ml</td><td>NM</td></tr><tr><td>OxiRed Probe (in DMSO)</td><td>0.2 ml</td><td>Red</td></tr><tr><td>HRP (lyophilized)</td><td>1 vial</td><td>Green</td></tr><tr><td>H<sub>2</sub>O<sub>2</sub> Substrate (0.88 M)</td><td>25 μl</td><td>Yellow</td></tr><tr><td>Stop Solution</td><td>1 ml</td><td>White</td></tr><tr><td>Catalase Positive Control</td><td>2 μl</td><td>Blue</td></tr></tbody></table>  | Components | Quantity | Color Code | Catalase Assay Buffer | 25 ml | NM | OxiRed Probe (in DMSO) | 0.2 ml | Red | HRP (lyophilized) | 1 vial | Green | H <sub>2</sub> O <sub>2</sub> Substrate (0.88 M) | 25 μl | Yellow | Stop Solution | 1 ml | White | Catalase Positive Control | 2 μl | Blue |
| Components                                       | Quantity  | Color Code |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Catalase Assay Buffer                            | 25 ml   | NM         |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| OxiRed Probe (in DMSO)                           | 0.2 ml  | Red        |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| HRP (lyophilized)                                | 1 vial  | Green      |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| H <sub>2</sub> O <sub>2</sub> Substrate (0.88 M) | 25 μl   | Yellow     |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Stop Solution                                    | 1 ml  | White      |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Catalase Positive Control                        | 2 μl  | Blue       |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Storage and Reagents Preparation                 | <p>Store the kit at -20°C, protect from light. Warm Assay Buffer to room temperature before use. Briefly centrifuge vials prior to opening. Read the entire protocol before performing the assay.</p> <ul style="list-style-type: none"><li>• <b>OxiRed Probe:</b> Briefly warm to completely melt the DMSO solution. Store at 4°C, protected from light. Use within two months.</li><li>• <b>HRP:</b> Dissolve with 220 μl Assay Buffer. Store at 4°C. Use within two months.</li><li>• <b>Positive Control Solution:</b> Add 500 μl Assay Buffer to Positive Control. Aliquot and store at -20°C. Diluted Positive Control solution is stable for 2-3 days at 4°C, for 2 months at -20°C.</li></ul> <p>Note: Keep samples, HRP and Catalase on ice while in use.</p>  |            |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |
| Assay Protocol                                   | <p><b>1. Sample and Positive Control Preparations:</b> Homogenize 0.1 gram tissues, or 10<sup>6</sup> Cells, or 0.2 ml Erythrocytes on ice in 0.2 ml cold Assay Buffer; Centrifuge at 10,000 x g for 15 minutes at 4°C; Collect the supernatant for assay, keep on ice. Liquid samples can be tested directly. Store samples at -80°C to assay later.</p> <p>Add 2 - 78 μl of samples or 1 - 5 μl Positive Control Solution into each well, and adjust volume to total 78 μl with Assay Buffer. Prepare sample High Control (HC) with the same amount of sample in separate wells then bring total volume to 78 μl with Assay Buffer. Add 10 μl of Stop Solution into the sample HC, mix and incubate at 25°C for 5 minutes to completely inhibit the catalase activity in samples as High Control. For unknown samples, we suggest testing several doses of your sample to ensure the readings are within the linear range.</p> <p>Reducing agents in samples interfere with the assay. Keep DTT or β-ME below 5 μM.</p> <p><b>2. H<sub>2</sub>O<sub>2</sub> Standard Curve:</b> Dilute 5 μl of 0.88M H<sub>2</sub>O<sub>2</sub> into 215 μl dH<sub>2</sub>O to generate 20 mM H<sub>2</sub>O<sub>2</sub>, then take 50 μl of the 20 mM H<sub>2</sub>O<sub>2</sub> and dilute into 0.95 ml dH<sub>2</sub>O to generate 1 mM H<sub>2</sub>O<sub>2</sub>. Add 0, 2, 4, 6, 8, 10 μl of 1 mM H<sub>2</sub>O<sub>2</sub> solution into 96-well plate to generate 0, 2, 4, 6, 8, 10 nmol/well H<sub>2</sub>O<sub>2</sub> standard. Bring the final volume to 90 μl with Assay Buffer. Add 10 μl Stop Solution into each well. For the fluorometric assay, dilute the standard H<sub>2</sub>O<sub>2</sub> 10-fold for the standard curve (0-1 nmol range).</p> <p>Note: Diluted H<sub>2</sub>O<sub>2</sub> is unstable, prepare fresh dilution each time.</p> <p><b>3. Catalase Reaction:</b> Add 12 μl fresh 1 mM H<sub>2</sub>O<sub>2</sub> into each well (samples, positive control, and sample HC) to start the reaction, incubate at 25°C for 30 minutes, and then add 10 μl Stop Solution into each sample well (Sample, Positive Control; do not add Stop Solution to the HC) to stop the reaction (Note: High Control and standard curve wells already contain Stop Solution).</p> <p><b>4. Develop Mix:</b> Mix enough reagents for the number of assays to be performed. For each well prepare a 50 μl Developer Mix containing:</p> |            |          |            |                       |       |    |                        |        |     |                   |        |       |  |       |        |               |      |       |                           |      |      |

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

Add 50 µl of the Developer Mix to each test samples, controls, and standards. Mix well and incubate at 25°C for 10 minutes. Measure OD 570 nm in a plate reader. Note: For low amounts of catalase, you can either increase the incubation time prior to adding the Stop Solution or use the fluorometric method. For the fluorometric method, decrease the 1 mM H<sub>2</sub>O<sub>2</sub> amount to 1.5 µl and OxiRed Probe to 0.3 µl in the reaction; compensate the volume with Assay Buffer.

**5. Calculation:** Signal change by catalase in sample is  $\Delta A = A_{HC} - A_{sample}$ . A<sub>HC</sub> is the reading of sample High Control, A<sub>sample</sub> is the reading of sample in 30 minutes. Plot the H<sub>2</sub>O<sub>2</sub> Standard Curve. Apply the  $\Delta A$  to the H<sub>2</sub>O<sub>2</sub> standard curve to get B nmol of H<sub>2</sub>O<sub>2</sub> decomposed by catalase in 30 minutes reaction. Catalase activity can be calculated:

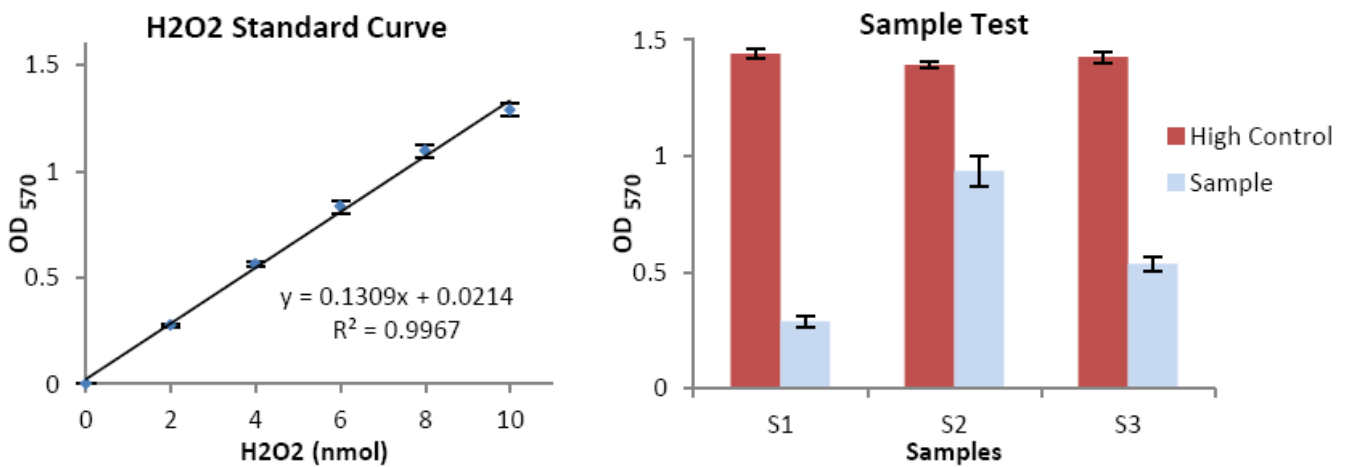
$$\text{Catalase Activity} = \frac{B}{30 \times V} \times \text{Sample Dilution Factor} = \text{nmol/min/ml} = \text{mU/mL}$$

Where: B is the decomposed H<sub>2</sub>O<sub>2</sub> amount from H<sub>2</sub>O<sub>2</sub> Standard Curve (in nmol).  
 V is the pretreated sample volume added into the reaction well (in ml).  
 30 is the reaction time 30 minutes.

Unit definition: One unit of catalase is the amount of catalase that decomposes 1.0 µmol of H<sub>2</sub>O<sub>2</sub> per minute at pH 4.5 at 25 °C.

Intended Use

For in vitro research use only. Not for diagnostic or therapeutic procedures.



**PromoCell GmbH**

Sickingenstr. 63/65  
 69126 Heidelberg  
 Germany

Email: info@promokine.info  
 www.promokine.info

**USA/Canada**

Phone: 1 – 866 – 251 – 2860 (toll free)  
 Fax: 1 – 866 – 827 – 9219 (toll free)

**Deutschland**

Telefon: 0800 – 776 66 23 (gebührenfrei)  
 Fax: 0800 – 100 83 06 (gebührenfrei)

**France**

Téléphone: 0800 90 93 32 (ligne verte)  
 Téléfax: 0800 90 27 36 (ligne verte)

**United Kingdom**

Phone: 0800 – 96 03 33 (toll free)  
 Fax: 0800 – 169 85 54 (toll free)

**Other Countries**

Phone: +49 6221 – 649 34 0  
 Fax: +49 6221 – 649 34 40