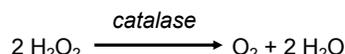


## EnzyChrom™ Catalase Assay Kit (ECAT-100)

### Quantitative Colorimetric/Fluorimetric Catalase Determination

#### DESCRIPTION

CATALASE (EC 1.11.1.6), is an ubiquitous antioxidant enzyme that catalyzes the decomposition of hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) to water and oxygen.



By preventing excessive H<sub>2</sub>O<sub>2</sub> build up, catalase allows important cellular processes which produce H<sub>2</sub>O<sub>2</sub> as a byproduct to occur safely. Deficiency in catalase activity has been associated with grey hair and peroxisomal disorder acatalasia. Simple, direct and high-throughput assays for catalase activity find wide applications. BioAssay Systems' improved assay directly measures catalase degradation of H<sub>2</sub>O<sub>2</sub> using a redox dye. The change in color intensity at 570nm or fluorescence intensity ( $\lambda_{em/ex} = 585/530\text{nm}$ ) is directly proportional to the catalase activity in the sample.

#### KEY FEATURES

**Sensitive and accurate.** Use 10  $\mu\text{L}$  sample. Linear detection range 0.2 to 5 U/L catalase activity.

**Simple and Convenient.** The procedure involves adding a Substrate to sample, incubation for 30 min, followed by a Detection Reagent and reading the optical density or fluorescence intensity.

#### APPLICATIONS:

**Direct Assays:** catalase activity in biological samples e.g. serum, plasma, urine, saliva, cell culture etc.

**Drug Discovery/Pharmacology:** effects of drugs on catalase activity.

#### KIT CONTENTS:

**Assay Buffer:** 25 mL    **HRP Enzyme:** 120  $\mu\text{L}$     **Dye Reagent:** 120  $\mu\text{L}$   
**H<sub>2</sub>O<sub>2</sub> Solution:** 100  $\mu\text{L}$  3% H<sub>2</sub>O<sub>2</sub>    **Positive Control:** 8  $\mu\text{L}$  Catalase

**Storage conditions.** The kit is shipped on ice. Store all components at -20°C. Shelf life of 6 months after receipt.

**Precautions:** reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### SAMPLE PREPARATION

Tissue (10 mg) and cells (10<sup>6</sup>) are homogenized in 200  $\mu\text{L}$  cold PBS. Centrifuge 10 min at 14,000 rpm to pellet any debris. Use clear supernatant for assay.

*Note: SH-containing reagents (e.g.  $\beta$ -mercaptoethanol, dithiothreitol) are known to interfere in this assay and should be kept below 10  $\mu\text{M}$  in the sample.*

#### ASSAY PROCEDURE

1. **Reagent Preparation.** Equilibrate all components to room temperature. Briefly centrifuge all tubes before opening. Keep thawed HRP Enzyme on ice.

For colorimetric assays, use a clear flat-bottom 96-well plate. For fluorimetric assays, use a solid black flat-bottom 96-well plate.

**Samples and Controls:** transfer 10  $\mu\text{L}$  sample into wells of the 96-well plate. In addition, for each assay run, prepare one sample blank well that contains only 10  $\mu\text{L}$  Assay Buffer.

Add 400  $\mu\text{L}$  Assay Buffer to Positive Control tube and mix well. Transfer 10  $\mu\text{L}$  of the reconstituted Positive Control into separate wells.

*Note: (1). For unknown samples, perform several dilutions to ensure that catalase activity is within the linear range 0.2 to 5 U/L. (2) The provided catalase serves as a positive control to ensure assay is working and should not be used to calculate the Sample catalase activity.*

2. **Enzyme Reaction.** Mix 5  $\mu\text{L}$  3% H<sub>2</sub>O<sub>2</sub> and 914  $\mu\text{L}$  dH<sub>2</sub>O (final 4.8 mM). Prepare enough 50  $\mu\text{M}$  H<sub>2</sub>O<sub>2</sub> Substrate for sample, positive control and sample blank by mixing, for each well, 1  $\mu\text{L}$  of the 4.8 mM H<sub>2</sub>O<sub>2</sub> with

95  $\mu\text{L}$  Assay Buffer. *Note: diluted H<sub>2</sub>O<sub>2</sub> is not stable. Prepare fresh dilutions for each experiment.*

Add 90  $\mu\text{L}$  of the 50  $\mu\text{M}$  Substrate to these wells to initiate the catalase reaction. Tap plate quick to mix. Incubate 30 min at room temperature. During the incubation time, proceed with Steps 3 and 4 below.

3. **H<sub>2</sub>O<sub>2</sub> Standard Curve.** Mix 40  $\mu\text{L}$  of the 4.8 mM H<sub>2</sub>O<sub>2</sub> with 440  $\mu\text{L}$  dH<sub>2</sub>O to yield 400  $\mu\text{M}$  H<sub>2</sub>O<sub>2</sub>. Prepare standards as shown in the Table below. Transfer 10  $\mu\text{L}$  standards into separate wells of the 96-well plate. Add 90  $\mu\text{L}$  Assay Buffer to the standards.

No	400 $\mu\text{M}$ H <sub>2</sub> O <sub>2</sub> + H <sub>2</sub> O	Vol ( $\mu\text{L}$ )	H <sub>2</sub> O <sub>2</sub> ( $\mu\text{M}$ )
1	100 $\mu\text{L}$ + 0 $\mu\text{L}$	100	400
2	60 $\mu\text{L}$ + 40 $\mu\text{L}$	100	240
3	30 $\mu\text{L}$ + 70 $\mu\text{L}$	100	120
4	0 $\mu\text{L}$ + 100 $\mu\text{L}$	100	0

4. **Detection.** Prepare enough Detection Reagent by mixing, for each reaction well (Sample, Control and Standard wells), 102  $\mu\text{L}$  Assay Buffer, 1  $\mu\text{L}$  Dye Reagent and 1  $\mu\text{L}$  HRP Enzyme.

At the end of the 30 min incubation (Step 2), add 100  $\mu\text{L}$  Detection Reagent per well. Tap plate to mix. Incubate for 10 min.

5. Read optical density at 570nm (550 to 585nm) or fluorescence intensity at  $\lambda_{em/ex} = 585/530\text{nm}$ .

#### CALCULATION

Subtract blank value (#4) from the standard values and plot the  $\Delta\text{OD}$  or  $\Delta\text{F}$  against standard concentrations. Determine the slope and calculate the catalase activity of Sample,

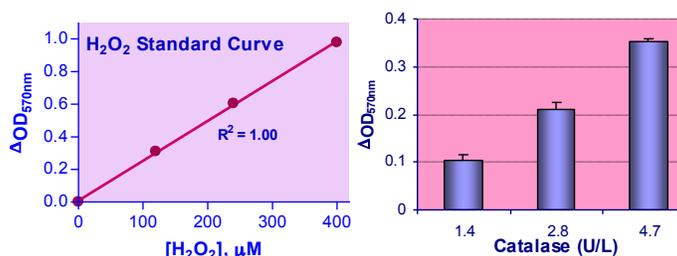
$$\text{Catalase (U/L)} = \frac{R_{\text{Sample Blank}} - R_{\text{Sample}}}{\text{Slope } (\mu\text{M}^{-1}) \times 30 \text{ min}} \times n$$

$R_{\text{SAMPLE Blank}}$  and  $R_{\text{SAMPLE}}$  are optical density or fluorescence intensity readings of the Sample Blank and Sample, respectively. Slope is determined from the standard curve. 30 min is the catalase reaction time.  $n$  is the sample dilution factor.

**Unit definition:** one unit is the amount of catalase that decomposes 1  $\mu\text{mole}$  of H<sub>2</sub>O<sub>2</sub> per min at pH 7.0 and room temperature.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, flat-bottom 96-well plates, plate reader.



#### PUBLICATIONS

- Chiu CC et al (2012). Beneficial Effects of Ocimum gratissimum Aqueous Extract on Rats with CCl<sub>4</sub>-Induced Acute Liver Injury. Evid Based Complement Alternat Med 2012:736752.
- Hadzi-Petrushev N et al (2012). L-2-oxothiazolidine-4-carboxylate influence on age- and heat exposure-dependent peroxidation in rat's liver and kidney. J Therm Biol 37(5):361-365.
- Zhu T et al (2012) Effects of the iron-chelating agent deferoxamine on triethylene glycol dimethacrylate, 2-hydroxyethyl methacrylate, hydrogen peroxide-induced cytotoxicity. J Biomed Mater Res B Appl Biomater 100(1):197-205.