

# Product Information & ELISA Manual

SOD (Superoxide Dismutase) Activity Assay Kit (Colorimetric) NBP3-24485

Enzyme-linked Immunosorbent Assay for quantitative detection.

#### Contact

Web: www.bio-techne.com/brands/novus-biologicals/ Email: nb-customerservice@bio-techne.com P: 888.506.6887 // P: 303.730.1950 // F: 303.730.1966 Novus kits are guaranteed for 6 months from date of receipt. For research use only. Not for diagnostic or therapeutic procedures.

# Table of contents

Assay summary	
Intended use	
Detection principle	4
Kit components & storage	5
Materials prepared by users	5
Reagent preparation	6
Sample preparation	7
The key points of the assay	9
Operating steps	
Calculation	
Appendix I Performance Characteristics	13
Appendix П Example Analysis	14
Statement	16

# Assay summary



# Intended use

This kit can be used to measure T-SOD, CuZn-SOD, Mn-SOD activity in serum, plasma, urine, cells, cell culture supernatant and tissue homogenate samples.

# **Detection principle**

Superoxide anion (O2•-) produced by xanthine and xanthine oxidase system can oxidize hydroxylamine to form nitrite which appear purplish red after chromogenic reaction. The SOD in the sample has a specific inhibitory effect on superoxide anion (O2•-), can reduce the content of nitrite. The OD value is lower than control and the activity of SOD is calculated through the formula. There are two kinds of SOD (CuZn-SOD, Mn-SOD) in the cells of higher animals, and the sum of them is equal to the total SOD. The activity of Mn-SOD in the sample will lost after sample pretreatment, but the activity of CuZn-SOD will not.



# Kit components & storage

Item	Component	Size 1 (50Assays)	Size 2 (100Assays)	Storage
Reagent 1	Buffer Solution	6 mL ×1 vial	12 mL ×1 vial	2-8°C, 12 months
Reagent 2	Nitrosogenic Agent	6 mL ×1 vial	12 mL ×1 vial	2-8°C, 12 months
Reagent 3	Substrate Solution	6 mL ×1 vial	12 mL ×1 vial	2-8°C, 12 months
Reagent 4	Enzyme Stock Solution	0.3 mL ×1 vial	0.6 mL ×1 vial	-20°C, 12 months
Reagent 5	Enzyme Diluent	6 mL ×1 vial	12 mL ×1 vial	2-8°C, 12 months
Reagent 6	Chromogenic Agent A	Powder ×1 vial	Powder ×1 vial	2-8°C, 12 months shading light
Reagent 7	Chromogenic Agent B	Powder ×1 vial	Powder ×1 vial	2-8°C, 12 months shading light
Reagent 8	Chromogenic Agent C	30 mL ×1 vial	60 mL ×1 vial	2-8°C, 12 months
Reagent 9	Extracting Solution	12.5 mL ×1 vial	25 mL ×1 vial	2-8°C, 12 months shading light

Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other. For a small volume of reagents, please centrifuge before use, so as not to obtain sufficient amount of reagents.

# Materials prepared by users

#### Instruments:

Spectrophotometer (550 nm), Micropipettor, Centrifuge, Incubator, Vortex mixer

#### **Reagents:**

Double distilled water, Normal saline (0.9% NaCl)

# **Reagent preparation**

- (1) Keep enzyme stock solution on ice during use. Equilibrate other reagents to room temperature before use.
- (2) The preparation of buffer working solution : For each tube, prepare 1 mL of buffer working solution (mix well 100 uL of buffer solution and 900 uL of double distilled water). Store at 2-8°C for 3 months.
- (3) The preparation of enzyme stock working solution : Before testing, please prepare sufficient enzyme stock working solution according to the test wells. For example, prepare 100 µL of enzyme stock working solution (mix well 5 uL of enzyme stock solution and 95 uL of enzyme diluent). Store at 2-8°C for 3 days.
- (4) The preparation of chromogenic agent A application solution : Dissolve one vial of chromogenic agent A with 70-80°C double distilled water to a final volume of 90 mL. Store at 2-8°C for 3 months protected from light.
- (5) The preparation of chromogenic agent B application solution : Dissolve one vial of chromogenic agent B with double distilled water to a final volume of 90 mL. Store at 2-8°C for 1 months protected from light.
- (6) The preparation of chromogenic agent : For each tube, prepare 2 mL of chromogenic agent (mix well 750 uL of chromogenic agent A application solution, 750 uL of chromogenic agent B application solution and 500 uL of chromogenic agent C). The chromogenic agent should be prepared on spot. Store at 2-8°C for 1 month protected from light.

# Sample preparation

# **1** Sample preparation

**Serum and plasma:** detect directly. If not detected on the same day, the serum or plasma can be stored at -80°C for a month.

**Urine:** Collect fresh urine and centrifuge at 10000 g for 15 minutes to remove insoluble material. Collect supernatant and keep it on ice for detection. If not detected on the same day, the urine can be stored at -80°C for a month.

#### **Tissue sample:**

- Harvest the amount of tissue needed for each assay (initial recommendation 20 mg).
- (2) Wash tissue in cold PBS (0.01 M, pH 7.4).
- (3) Homogenize 20 mg tissue in 180 µL normal saline (0.9% NaCl) with a dounce homogenizer at 4°C.
- (4) Centrifuge at 10000×g for 10 min at 4°C to remove insoluble material. Collect supernatant and keep it on ice for detection.
- Meanwhile, determine the protein concentration of supernatant (E-BC-K318-M).

#### Cells:

- Harvest the number of cells needed for each assay (initial recommendation 10<sup>6</sup> cells).
- (2) Wash cells with PBS (0.01 M, pH 7.4).
- (3) Homogenize 10<sup>6</sup> cells in 300-500 μL normal saline (0.9% NaCl) with a ultrasonic cell disruptor at 4°C.
- (4) Centrifuge at 10000×g for 10 min at 4°C to remove insoluble material. Collect supernatant and keep it on ice for detection.
- Meanwhile, determine the protein concentration of supernatant (E-BC-K318-M).

## **②** Dilution of sample

The optimal sampling volume are different for different species, the SOD also are different for different samples. It is recommended to take  $2\sim3$  samples to do a pre-experiment, diluting a series of diluent and determine the dilution factor when the SOD inhibition ratio is  $15\%\sim55\%$  (the optimal inhibition ratio is the range of  $25\%\sim45\%$ .) before formal experiment.

The recommended dilution factor for different samples is as follows (for reference only):

Sample type	Dilution factor	The volume added to the
		reaction (µL)
Mouse serum	3-5	20-30
Human serum	1	30-40
Rat serum	3-5	20-30
Rat plasma	1-2	20-30
Human hydrothorax	1	30-50
Human urine	1	30-50
10% Mouse liver tissue homogenate	50-80	20-30
10% Mouse brain tissue homogenate	8-12	20-30
10% Mouse kidney tissue homogenate	10-20	20-30
HepG2 cells (3.27 mgprot/mL)	5-10	20-30

Note: The diluent is normal saline (0.9% NaCl). For the dilution of other sample types, please do pretest to confirm the dilution factor

## The key points of the assay

- Determine optimal sampling volume of each sample before formal experiment. Calculate the inhibition ratio of serial sampling volume, and choose the optimal sampling volume when inhibition ratio in the range of 25%~45%.
- 2 The optimal sampling volume are different for different species, the SOD also are different for different samples. So it is best to do a pre-test to determining optimal sampling volume for a new sample.
- (3) It is best to reserve 3 paralleled tubes with different sampling volumes in pretest for determining the optimal sampling volume.
- (4) Adjust sampling volume: If inhibition ratio >55%, need to dilute the sample or decrease the sampling volume than take the test. If inhibition ratio < 15%, need to increase the sampling volume.</p>
- (5) There should be no bubbles in the wells of the microplate when measuring the OD value.
- (6) The multichannel pipeter is recommended to shorten the time and reduce the error between wells.

# **Operating steps**

#### Sample pretreatment

- Take 0.1 mL sample and add 0.1 mL extracting solution. Mix thoroughly with a vortex mixer for 1 min by vortex mixer and centrifuge at 3500 g for 15 min. Take supernatant for CuZn-SOD measurement.
- (2) Take 0.1 mL normal saline and add 0.1 mL extracting solution. Mix thoroughly for 1 min with a vortex mixer and centrifuge at 3500 g for 15 min. Take supernatant as control of CuZn-SOD.

#### The measurement of samples

 Sample tube of T-SOD: take 1 mL of buffer working solution and a\* μL of sample to the corresponding tubes.

Control tube of T-SOD: take 1 mL of buffer working solution and  $a^* \mu L$  of double distilled water to the corresponding tubes.

Sample tube of CuZn-SOD: take 1 mL of buffer working solution and  $a^* \mu L$  of supernatant for CuZn-SOD to the corresponding tubes.

Control tube of CuZn-SOD: take 1 mL of buffer working solution and  $a^* \mu L$  of supernatant as control of CuZn-SOD to the corresponding tubes.

- (2) Add 0.1 mL of nitrosogenic agent, 0.1 mL of substrate solution, 0.1 mL of enzyme stock working solution to each tube.
- (3) Mix fully with a vortex mixer and incubate at 37°C for 40 min. (When the room temperature is below 20°C, the time of incubation can be extended to 45 min.).
- (4) Add 2 mL of chromogenic agent to each tube.
- (5) Mix fully and stand for 10 min at room temperature. Set the spectrophotometer to zero with double distilled water and measure the OD values of each tube at 550 nm with 1 cm optical path cuvette.

# Calculation

# The sample:

# 1. Serum (plasma), cell culture medium and other liquid samples:

**Definition:** When SOD inhibition ratio in 1 mL reaction system reach 50%, the corresponding enzyme level is 1 SOD activity unit (U).

$$\frac{\text{T-SOD activity}}{(U/mL)} = i_1 \div 50\% \times \frac{V_1}{V_2} \times f$$

$$\frac{\text{CuZn-SOD activity}}{(U/mL)} = i_2 \div 50\% \times \frac{V_1}{V_2} \times f$$

# 2. Tissue and cell samples:

**Definition:** When SOD inhibition ratio of 1 mg of tissue protein in 1 mL reaction system reach 50%, the corresponding enzyme level is 1 SOD activity unit (U)

$$\begin{array}{rl} \text{T-SOD activity}\\ (U/\text{mgprot}) &= i_1 \stackrel{\div}{\to} 50\% \times \frac{V_1}{V_2} \times f \stackrel{\div}{\to} C\\ \text{CuZn-SOD activity}\\ ((U/\text{mgprot}) &= i_2 \stackrel{\div}{\to} 50\% \times \frac{V_1}{V_2} \times f \stackrel{\div}{\to} C_{\text{pr}} \end{array}$$

Mn-SOD activity =T-SOD activity -CuZn-SOD activity

# [Note]

i1: Inhibition ratio of T-SOD.

i2: Inhibition ratio of CuZn-SOD.

 $A_1$ : The OD value of T-SOD<sub>Control</sub>.

A<sub>2</sub>: The OD value of T-SOD<sub>Sample</sub>.

A<sub>3</sub>: The OD value of CuZn-SOD<sub>Control</sub>.

A4: The OD value of CuZn-SOD<sub>Sample</sub>.

V<sub>1</sub>: The total volume of the reaction system (mL).

V<sub>2</sub>: The volume of sample added to the reaction system (mL).

f: Dilution factor of sample before test..

C<sub>pr</sub>: Concentration of protein in sample, mgprot/mL.

# **Appendix I Performance Characteristics**

#### 1. Parameter:

#### **Intra-assay Precision**

Three human serum samples were assayed in replicates of 20 to determine precision within an assay (CV = Coefficient of Variation).

Parameters	Sample 1	Sample 2	Sample 3
Mean (U/mL)	6.70	58.50	132.50
%CV	4.3	3.9	3.8

#### **Inter-assay Precision**

Three human serum samples were assayed 20 times in duplicate by three operators to determine precision between assays.

Parameters	Sample 1	Sample 2	Sample 3
Mean (U/mL)	6.70	58.50	132.50
%CV	7.1	6.9	7.6

## Recovery

Take three samples of high concentration, middle concentration and low concentration to test the samples of each concentration for 6 times parallelly to get the average recovery rate of 96%.

	Sample 1	Sample 2	Sample 3
Expected Conc. (U/mL)	18.5	67.5	123
Observed Conc. (U/mL)	18.3	63.5	116.9
Recovery rate (%)	99	94	95

#### Sensitivity

The analytical sensitivity of the assay is 2.03 U/mL. This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 20 times, and calculating the corresponding concentration.

### **Appendix II Example Analysis**

#### **Example analysis :**

**The detection of T-SOD:** Take 10% mouse heart tissue homogenate, dilute for 10 times with normal saline (0.9% NaCl), then take 20  $\mu$ L of diluted sample, and carry the assay according to the operation steps. The results are as follows:

The average OD value of T-SOD<sub>Sample</sub> is 0.252, the average OD value of T-SOD<sub>Control</sub> is 0.522, the concentration of protein in 10% mouse heart tissue homogenate is 6.84 mgprot/mL, and the calculation result is:

$$\frac{\text{T-SOD activity}}{(\text{U/mgprot})} = \frac{0.522 \cdot 0.252}{0.522} \div 50\% \times \frac{3.02}{0.02} \times 10 \div 6.84 = 228.37 \text{ U/mgprot}$$

**The detection of CuZn-SOD:** pretreat the diluted sample with extracting solution, and carry the assay according to the operation steps. The results are as follows: The average OD value of CuZn-SOD<sub>Sample</sub> is 0.365, the average OD value of CuZn-SOD<sub>Control</sub> is 0.571, the concentration of protein in 10% mouse heart tissue homogenate is 6.84 mgprot/mL, and the calculation result is:

$$\frac{\text{CuZn-SOD activity}}{(\text{U/mgprot})} = \frac{0.571 - 0.365}{(\frac{0.571}{0.571}) \div 50\% \times \frac{3.02}{0.02} \times 10 \div 6.84 = 159.29 \text{ U/mgprot}}$$

Detect human serum (V<sub>2</sub>=30  $\mu$ L), mouse serum (dilute for 4 times, V<sub>2</sub>=20  $\mu$ L), 10% mouse heart tissue homogenate (the concentration of protein is 6.84 mgprot/mL, dilute for 10 times, V<sub>2</sub>=20  $\mu$ L) and HepG2 cells (the concentration of protein is 3.27 mgprot/mL, dilute for 10 times, V<sub>2</sub>=25  $\mu$ L) according to the protocol, the result is as follows :





#### Statement

- 1. This assay kit is for Research Use Only. We will not response for any arising problems or legal responsibilities causing by using the kit for clinical diagnosis or other purpose.
- 2. Please read the instructions carefully and adjust the instruments before the experiments. Please follow the instructions strictly during the experiments.
- 3. Protection methods must be taken by wearing lab coat and latex gloves.
- 4. If the concentration of substance is not within the detection range exactly, an extra dilution or concentration should be taken for the sample.
- 5. It is recommended to take a pre-test if your sample is not listed in the instruction book.
- 6. The experimental results are closely related to the situation of reagents, operations, environment and so on. Novus will guarantee the quality of the kits only, and NOT be responsible for the sample consumption caused by using the assay kits. It is better to calculate the possible usage of sample and reserve sufficient samples before use.