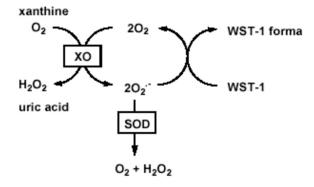
#### Technical Bulletin

# Superoxide Dismutase (SOD) Activity Kit

#### Catalog Number MAK379

# **Product Description**

Superoxide dismutase (SOD) is one of the most important antioxidative enzymes. It catalyzes the dismutation of the superoxide anion into hydrogen peroxide and molecular oxygen. The sensitive SOD assay kit utilizes WST-1 that produces a water-soluble formazan dye upon reduction with superoxide anion. The rate of the reduction with a superoxide anion is linearly related to the xanthine oxidase (XO) activity and is inhibited by SOD (below). Therefore, the inhibition activity of SOD can be determined by a colorimetric method.



The kit is suitable for the determination of superoxide dismutase (SOD) activity in blood, tissue and cell lysates.

# **Components**

The kit is sufficient for 100 colorimetric assays in 96-well plates.

•	WST Solution Catalog Number MAK379A	1 mL
•	SOD Enzyme Solution Catalog Number MAK379B	20 μL
•	SOD Assay Buffer Catalog Number MAK379C	20 mL
•	SOD Dilution Buffer Catalog Number MAK379D	10 mL

# Reagents and Equipment Required but Not Provided

- Pipetting devices and accessories (including multichannel pipettor)
- 96-well flat-bottom plate. It is recommended to use clear plates for colorimetric assays. Cell culture or tissue culture treated plates are **not** recommended.
- Spectrophotometric multiwell plate reader
- Refrigerated microcentrifuge capable of RCF  $\geq$ 14,000  $\times$  g
- Dounce tissue grinder set (Catalog Number D9063 or equivalent)



- Mitochondria/Cytosol Fractionation Kit (Catalog Number MIT1000)
- Phosphate Buffered Saline (Catalog Number P3813 or equivalent)
- Trizma<sup>®</sup> hydrochloride (Catalog Number T3253 or equivalent)
- Triton™ X-100 (Catalog Number X100)
- 2-Mercaptoethanol (Catalog Number M6250 or equivalent)
- Phenylmethanesulfonyl fluoride (PMSF) (Catalog Number P7626)

#### **Precautions and Disclaimer**

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

# Storage/Stability

The kit is shipped on wet ice. Store components at 2-8 °C. Briefly centrifuge small vials at low speed prior to opening.

# **Preparation Instructions**

<u>WST Working Solution</u>: Dilute the 1 mL of WST solution with 19 mL of SOD Assay Buffer. The diluted solution is stable for up to 2 months at 2-8 °C.

Enzyme Working Solution: Centrifuge the SOD Enzyme Solution for 5 seconds. Mix well by pipetting (The step is necessary, as the enzyme has two layers and must be mixed well before dilution). Dilute 15  $\mu$ L of the mixed enzyme with 2.5 mL of SOD Dilution Buffer. The diluted enzyme solution is stable for up to 3 weeks at 2-8 °C.

#### Sample Preparation

Blood samples: Collect blood using citrate or EDTA. Centrifuge at  $1,000 \times g$  for 10 minutes at 2-8 °C. Transfer the plasma layer to a new tube without disturbing the buffy layer and store at -80 °C until ready for analysis. Remove the buffy layer from the red cell pellet. Resuspend the erythrocytes in  $5\times$  volume of ice-cold purified water and centrifuge at  $10,000\times g$  for 10 minutes to pellet the erythrocyte membranes. Store the supernatant at -80 °C until ready for analysis. Plasma can be diluted approximately  $3-10\times$  and the red cell lysate diluted approximately  $100\times$  prior to SOD assay.

<u>Tissue and cells</u>: Tissue should be perfused with PBS to remove any red blood cells. Homogenize tissue or lyse cells in ice-cold 0.1 M Tris-HCl, pH 7.4 containing 0.5% Triton X-100, 5 mM 2-mercaptoethanol, 0.1 mg/ml PMSF. Centrifuge the crude tissue homogenate/cell lysate at  $14000 \times g$  for 5 minutes at 2-8 °C and discard the cell debris. The supernatant contains total SOD activity from cytosolic and mitochondria.

#### **Procedure**

<u>Note</u>: If it is desired to measure SOD activity from cytosol and mitochondria separately, cytosol and mitochondria can be separated by using the Mitochondria/Cytosol Fractionation Kit. SOD activity is then measured from the Mitochondria and Cytosol fractions separately.

#### Assay Procedure

1. Refer to Table 1 for the preparation of sample(s) and blank reaction mixes. Note: A SOD standard is **not** included with the kit. If desired, set up wells for standards in the same manner as the sample. Since the superoxide will release immediately after the addition of Enzyme Working Solution to each well, use a multichannel pipettor to avoid reaction time lag of each well.



Table 1. Reaction Mix Preparation

Reagent	Sample (s)	Blank 1	Blank 2	Blank 3
Sample(s)	20 μL		20 μL	
Purified Water		20 μL		20 μL
WST Working Soln.	200 μL	200 μL	200 μL	200 μL
Enzyme Working Soln.	20 μL	20 μL		
Dilution Buffer			20 μL	20 μL

2. Mix well and incubate plates at 37 °C for 20 minutes.

# <u>Measurement</u>

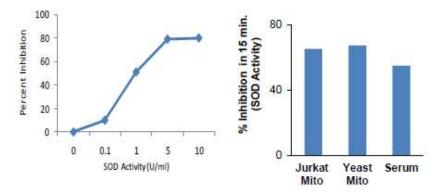
- 1. Read the absorbance at 450 nm using a microplate reader.
- 2. Calculate the SOD activity (inhibition rate %):

# **Results**

SOD Activity (inhibition rate %) =

$$(A_{450} \text{ blank1} - A_{450} \text{ blank3}) - (A_{450} \text{ sample} - A_{450} \text{ blank2}) \times 100$$
  
 $(A_{450} \text{ blank1} - A_{450} \text{ blank3})$ 

Figure 1.



SOD Activity (% inhibition rate): human serum (10  $\mu$ L) and isolated mitochondria from Jurkat cells (10  $\mu$ g), and yeast (*Saccromyces cerevisiae*, 100  $\mu$ g), was used to determine SOD Activity according to the kit protocol. Activity was measured at 15 minutes at 37 °C.



#### **Notice**

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

#### **Contact Information**

For the location of the office nearest you, go to SigmaAldrich.com/offices.

#### **Technical Service**

Visit the tech service page on our web site at SigmaAldrich.com/techservice.

#### **Standard Warranty**

The applicable warranty for the products listed in this publication may be found at <a href="SigmaAldrich.com/terms">SigmaAldrich.com/terms</a>.

MAK379 Technical Bulletin Rev 05/2021



The life science business of Merck KGaA, Darmstadt, Germany

operates as MilliporeSigma in the U.S. and Canada.

