

# SOD activity kit

ELS offers a large collection of superoxide dismutase products for oxidative stress research.

This is a versatile assay for the determination of SOD activity. Differentiate between the activity of three SOD isoenzymes: MnSOD, FeSOD, and Cu/ZnSOD. Obtain higher throughput sample analysis with a convenient colorimetric 96 well plate format and fast results in just 10 minutes.

SOD (Superoxide dismutase) is responsible for the elimination of cytotoxic active oxygen by catalyzing the dismutation of the superoxide radical to oxygen and hydrogen peroxide. There are three SOD isoenzymes in mammalian cells, they are: EC SOD (extracellular SOD), Cu/Zn SOD (copper and zinc-containing SOD) and Mn SOD (manganese-containing SOD).

Citations: 49

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-900-157

5x96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Versatile - differentiate between activity of three SOD isoenzymes: MnSOD, FeSOD, and Cu/ZnSOD
- Higher throughput - colorimetric 96 well plate format
- Fast - results in just 10 minutes

## Handling & Storage

**Use/Stability** Store all components at 4°C, except SOD standard and Xanthine Oxidase at -20°C.

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Superoxide dismutase

**Application** Activity assay, Colorimetric detection

**Application Notes** For the measurement of total SOD activity in cell lysates and tissue from any species. Cited sample type includes serum and gastric mucosa homogenates (ref.).

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Microtiter plate, Standard, SOD buffer concentrate, Xanthine oxidase, Xanthine solution, Triton x-100, WST-1 reagent



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)