

Superoxide Dismutase Assay

A plate based colorimetric assay to measure superoxide dismutase activity in various samples.

Superoxide is formed as a side product of a number of redox enzymes, such as mitochondrial complex I and complex II during respiration and NADPH oxidase as a part of the immune system. Superoxide is toxic due to its high reactivity and the level of superoxide must be controlled in living organisms. Superoxide dismutase (SOD) is an important antioxidative enzyme that catalyzes the dismutation, or removal, of superoxide into hydrogen peroxide and molecular oxygen. SODs are important because they are an organism's first defense against harmful superoxide anions. AkrivisBio's SOD Activity Assay provides a simple, quick and quantitative method to measure SOD activity in samples. In this assay, superoxide produced by the action of xanthine oxidase reduces WST from a nearly colorless tetrazolium to a highly colored formazan, which can be detected by absorbance at 450 nm. Superoxide is scavenged in an amount proportional to the SOD activity present in samples. Thus, greater the activity of SOD in samples, lesser is the formation of the colored formazan dye.

Ordering Information

[Order Online »](#)

AKR-MA-0108	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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