

Catalase Activity Assay

A plate based colorimetric assay to measure catalase activity in a wide range of biological samples.

Hydrogen peroxide is a common metabolic byproduct of many cellular processes. A potent oxidant, it's toxic due to its ability to damage cell structures, proteins and DNA. Catalase evolved to remove hydrogen peroxide and is one of the most efficient enzymes with turnover numbers approaching $3 \times 10^6 \text{ s}^{-1}$. It acts in maintaining a level of ROS in cells necessary for normal physiological functions without allowing harm from excessive amounts. AkrivisBio's Catalase Activity Assay is a simple, sensitive way to measure catalase activity in a range of biological samples. In the assay, any hydrogen peroxide not consumed by catalase is utilized by peroxidase to convert ADHP to resorufin with strong absorbance and fluorescence.

Ordering Information

[Order Online »](#)

AKR-MA-0143	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