

Catalase fluorometric detection kit

The Catalase fluorometric detection kit provides a higher throughput sensitive method for measuring catalase activity in just one hour. This is a homogeneous mix-and-read assay that enables you to monitor multiple time points for a kinetic read. Positive control is provided.

Catalase is an antioxidant enzyme that catalyses the decomposition of hydrogen peroxide (H_2O_2) to water and oxygen. Catalase is ubiquitously expressed in mammalian and non-mammalian aerobic cells containing the cytochrome system and functions to alleviate oxidative damage by catalyzing metabolism of the peroxide substrate. Accumulation of H_2O_2 can result in cellular damage through oxidation of proteins, DNA, and lipids, and has been linked to numerous pathologies including inflammation, cancer, diabetes, Parkinson's disease, and Alzheimer's.

Citations: 12

[View Online »](#)

Ordering Information

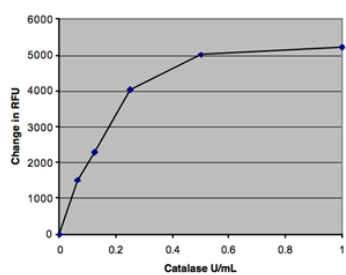
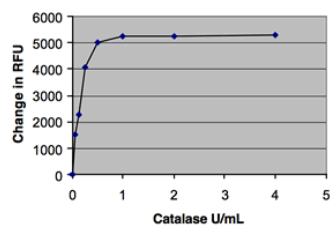
[Order Online »](#)

ADI-907-027	500 tests
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Homogeneous - mix-and-read format easily adaptable to HTS applications
- Sensitive - fluorescent detection system
- Fast - results in just 1 hour



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection, HTS

Application Notes For the measurement of catalase in cell lysates from any species.

Contents Reaction buffer concentrate, Detection Reagent, Horseradish peroxidase concentrate, Hydrogen peroxide, Catalase enzyme

Sensitivity (range 0- 4 U/ml)

Species Reactivity Species independent

UniProt ID P00432



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)