

Hydrogen Peroxide Assay

A plate based colorimetric/fluorometric assay to measure hydrogen peroxide in a wide range of samples.

Hydrogen peroxide is important in cellular metabolism due to its role as a reactive oxygen species. Although considered a harmful byproduct of metabolism, low levels of H₂O₂ are essential for some physiological processes. H₂O₂ acts as a signaling molecule, participating in signaling pathways that regulate responses to stress, growth, and immune functions. H₂O₂ is produced during aerobic respiration as a byproduct of cellular energy production. Mitochondria are a major source of H₂O₂ production. The controlled generation of H₂O₂ serves as an intermediary in cellular signaling, assisting in the activation of various enzymes involved in key metabolic pathways. H₂O₂ plays a role in the immune response, acting as a defense mechanism against invading pathogens. Immune cells use hydrogen peroxide to neutralize bacteria and viruses, highlighting its role in combating infections. AkrisBio's Hydrogen Peroxide Assay is a simple, sensitive means of measuring peroxide in a variety of samples. It can be used in either absorbance or fluorescence-based measurements.

Ordering Information

[Order Online »](#)

AKR-MA-0145	100 wells
--------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°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