

Hydrogen peroxide chemiluminescent detection kit

The Hydrogen Peroxide chemiluminescent detection kit provides a fast and higher throughput 96-well plate based homogeneous assay for the measurement of hydrogen peroxide. This kit offers more sensitivity over the Hydrogen Peroxide colorimetric detection kit (Cat. #ADI-907-015) but is not recommended for samples that contain phenol red or hemoglobin.

Hydrogen Peroxide (H_2O_2) is a reactive oxygen metabolic by-product that serves as a key regulator for a number of oxidative stress related states. Functioning through NFkB and other factors, Hydrogen peroxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis, diabetic vasculopathy, osteoporosis, and a number of neurodegenerative diseases.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

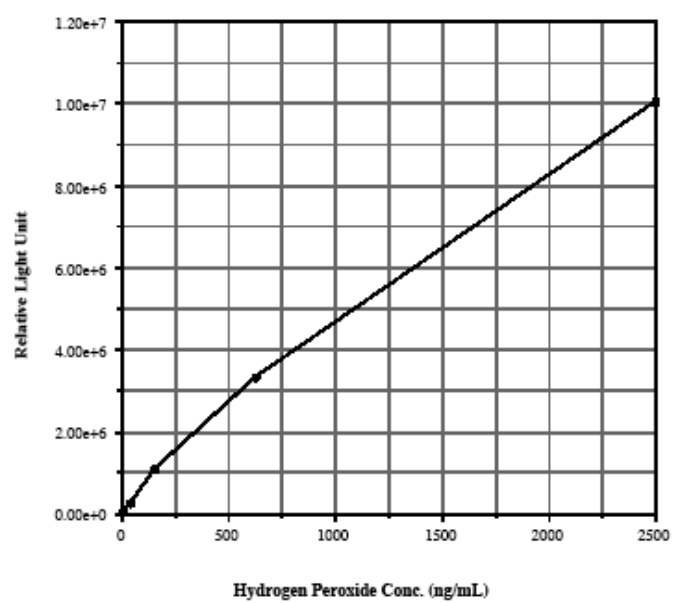
ADI-907-012	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Immediate results in < 10 minutes
- Extremely sensitive quantitation of hydrogen peroxide

Standard Curve in Buffer



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Chemiluminescence

Application Notes For the quantitative determination of hydrogen peroxide in culture supernatants from any species.

Contents White Microtiter plate, Standard, Substrate, Trigger solution

Sensitivity 2.39 ng/ml (range 2.44 – 2,500 ng/ml)

Species Reactivity Species independent



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)