

Hydrogen peroxide colorimetric detection kit

The Hydrogen Peroxide colorimetric detection kit provides a fast and higher throughput 96-well plate based homogeneous assay for the measurement of hydrogen peroxide in a wide variety of samples.

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: 26

[View Online »](#)

Ordering Information

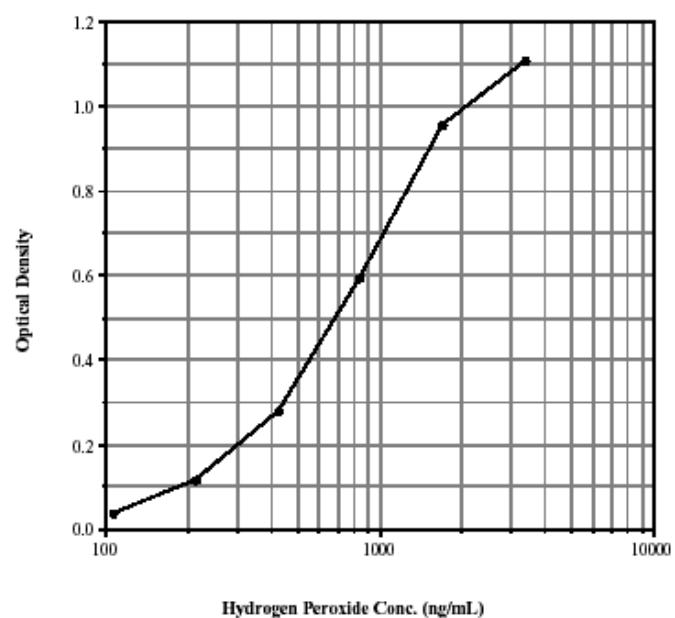
[Order Online »](#)

| | |
|-------------|----------|
| ADI-907-015 | 96 wells |
|-------------|----------|

Manuals, SDS & CofA

[View Online »](#)

- Fast results in just 30 minutes
- Sensitive to 1.51µM



Handling & Storage

Use/Stability Store all components at 4°C, except Color Reagent at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, HTS

Application Notes For the quantitative determination of hydrogen peroxide in culture supernatants, serum, plasma, and urine from any species. Cited sample type includes tissue (ref. 1).

Contents Half-Area Microtiter plate, Standard, Color reagent

Sensitivity 51.25 ng/ml (range 106.25 – 3,400 ng/ml)

Species Reactivity Species independent

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)