

Hydrogen peroxide fluorometric detection kit

The Hydrogen Peroxide fluorometric detection kit provides fast results in just 10 minutes for the measurement of either hydrogen peroxide or peroxidase activity in a convenient one-step homogeneous assay.

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 »](#)

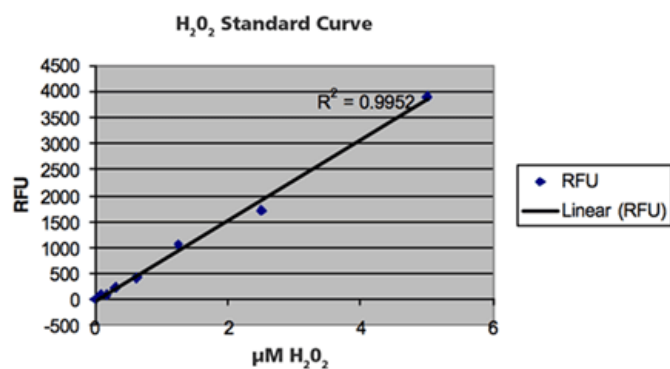
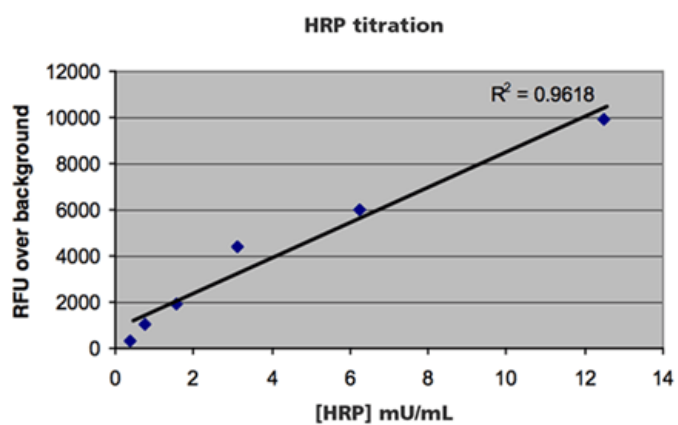
ADI-907-028

500 tests

Manuals, SDS & CofA

[View Online »](#)

- Flexible - measure hydrogen peroxide or peroxidase activity
- Higher throughput - one step homogeneous mix-and-read assay
- Fast - results in just 10 minutes



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 hydrogen peroxide in cell lysates and culture supernatants from any species.

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

Sensitivity (range 0 – 10 µM (H₂O₂); 0 – 25 mU/ml (peroxidase))

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)