## 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 NF $\kappa$ B 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** »

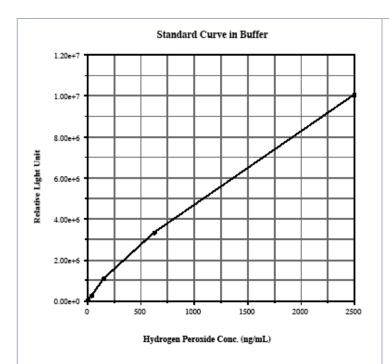
ADI-907-012

96 wells

Manuals, SDS & CofA

View Online »

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



## **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

Last modified: May 29, 2024



eu@enzolifesciences.com