## 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 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: 25

View Online »

**Ordering Information** 

**Order Online** »

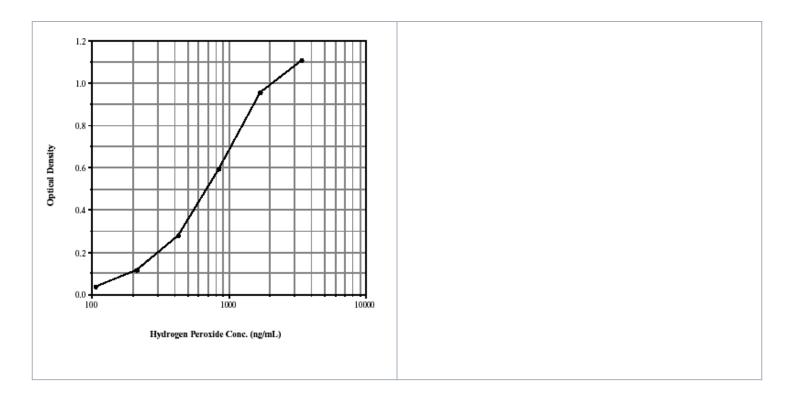
ADI-907-015

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

