

# Peroxide Assay Kit ab272537

[1 Image](#)

### Overview

|                         |   |
|-------------------------|---|
| <b>Product name</b>     | Peroxide Assay Kit  |
| <b>Detection method</b> | Colorimetric  |
| <b>Sample type</b>      | Urine, Serum, Plasma, Other biological fluids, Cell Lysate, Cell culture media  |
| <b>Assay type</b>       | Quantitative  |
| <b>Range</b>            | 0.2 $\mu$ M - 30 $\mu$ M  |
| <b>Product overview</b> | <p>Peroxide Assay Kit (ab272537) is a simple, direct and automation-ready procedures for quantitative determination of peroxide find wide applications in research and drug discovery. This assay kit is designed to measure peroxide concentration in biological samples without any pretreatment. The improved method utilizes the chromogenic <math>\text{Fe}^{3+}</math>-xylenol orange reaction, in which a purple complex is formed when <math>\text{Fe}^{2+}</math> provided in the reagent is oxidized to <math>\text{Fe}^{3+}</math> by peroxides present in the sample. The intensity of the color, measured at 540-610nm, is an accurate measure of the peroxide level in the sample. The optimized formulation substantially reduces interference by substances in the raw samples.</p> |

**Sensitive and accurate.** Enhanced color intensity using sorbitol. Detection range 0.2  $\mu$ M (7 ng/mL) to 30  $\mu$ M (1,020 ng/mL)  $\text{H}_2\text{O}_2$  in 96-well plate assay.

**Simple and high-throughput.** The procedure involves addition of a single detection reagent and incubation for 30 min. Can be readily automated as a high-throughput assay in 96-well plates for thousands of samples per day.

|                            |   |
|----------------------------|---|
| <b>Tested applications</b> | <b>Suitable for:</b> Functional Studies |
| <b>Platform</b>            | Microplate                              |

### Properties

|                             |   |
|-----------------------------|---|
| <b>Storage instructions</b> | Store at +4°C. Please refer to protocols. |
|-----------------------------|---|

| Components  | 250 tests |
|---|-----------|
| Reagent A   | 1 x 1ml   |
| Reagent B   | 1 x 50ml  |
| Standard (3% stabilized H <sub>2</sub> O <sub>2</sub> ) | 1 x 100µl |

## Relevance

Hydrogen Peroxide (H<sub>2</sub>O<sub>2</sub>) is a reactive oxygen metabolic byproduct that serves as a key regulator for a number of oxidative stress-related states. Functioning through NF-κB and other factors, hydroperoxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis, diabetic vasculopathy, osteoporosis, neuro-degenerative diseases, Down's syndrome and immune system diseases.

## Applications

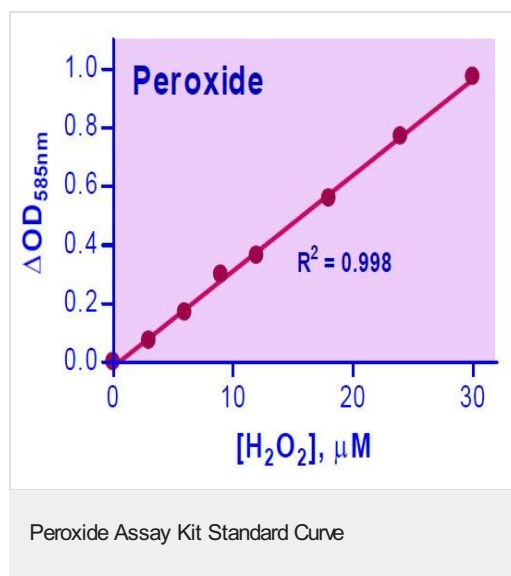
### The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab272537 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

| Application        | Abreviews | Notes                                    |
|--------------------|-----------|--|
| Functional Studies |           | Use at an assay dependent concentration. |

## Images



**Typical standard curve** – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet

- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors