

### Peroxide Assay Kit ab272537

画像数 1

医薬用外劇物

#### 製品の概要

製品名	Peroxide Assay Kit
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Plasma, Other biological fluids, Cell Lysate, Cell culture media
アッセイタイプ	Quantitative
検出範囲	0.2 $\mu$ M - 30 $\mu$ M
製品の概要	<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> <p><b>Sensitive and accurate.</b> Enhanced color intensity using sorbitol. Detection range 0.2 <math>\mu</math>M (7 ng/mL) to 30 <math>\mu</math>M (1,020 ng/mL) <math>\text{H}_2\text{O}_2</math> in 96-well plate assay.</p> <p><b>Simple and high-throughput.</b> 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.</p>
アプリケーション	<b>適用あり:</b> Functional Studies
試験プラットフォーム	Microplate

#### 製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	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

## 関連性

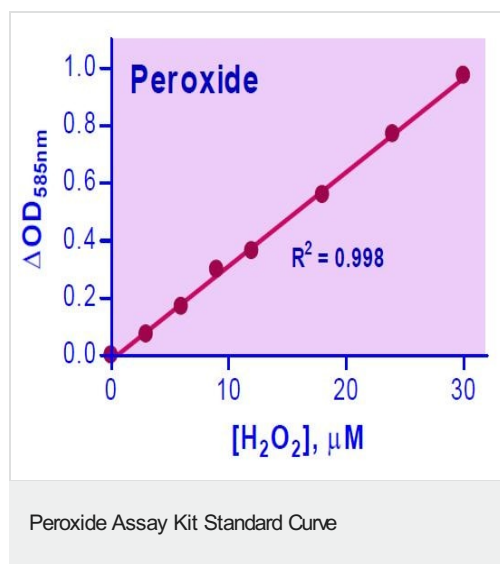
Hydrogen Peroxide (H<sub>2</sub>O<sub>2</sub>) is a reactive oxygen metabolic byproduct that serves as a key regulator for a number of oxidative stress-related states. Functioning through NF-κB and other factors, hydroperoxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis,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

## アプリケーション

**The Abpromise guarantee** Abpromise保証は、次のテスト済みアプリケーションにおけるab272537の使用に適用されます  
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

## 画像



**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.co.jp/abpromise> or contact our technical team.

#### **Terms and conditions**

---

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