

Hydrogen Peroxide Assay Kit (Cell-based) ab138874

★★★★★ [1 Abreviews](#) [5 References](#) [画像数 1](#)

製品の概要

製品名	Hydrogen Peroxide Assay Kit (Cell-based)
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based
検出感度	3 μ M
全工程の試験時間	0h 30m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Hydrogen Peroxide Assay Kit (Cell-based) (ab138874) uses our AbGreen Indicator to quantify hydrogen peroxide in live cells. The indicator is cell-permeable, and generates green fluorescence when it reacts with hydrogen peroxide.</p> <p>The assay is in a “mix and read” assay format that is compatible with HTS liquid handling instruments and a 96-well or 384-well plate format. It provides a sensitive, one-step fluorometric assay to detect as little as 0.3 nanomoles of H_2O_2 in a 100 μL assay volume (3 μM). Results can be read by either a fluorescence microplate reader at Ex/Em = 490/520 nm for H_2O_2 detection in solution or a fluorescence microscopy for live cell H_2O_2 imaging.</p> <p>Hydrogen peroxide assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add green indicator solution - analyze with microplate reader for 15-60 min. <p>Alternatively, for live cell assay:</p> <ul style="list-style-type: none"> - add green indicator solution to cells - incubate for 5-60 min whilst analyzing with microplate reader or fluorescence microscope
特記事項	<p>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. It is involved in many biological events that are linked to asthma, atherosclerosis, diabetic vasculopathy, osteoporosis, a number of neurodegenerative diseases and Down’s syndrome. The measurement of this reactive species is helpful for determining how oxidative stress modulates various intracellular pathways.</p>
試験プラットフォーム	Fluorescence microscope

製品の特性

保存方法

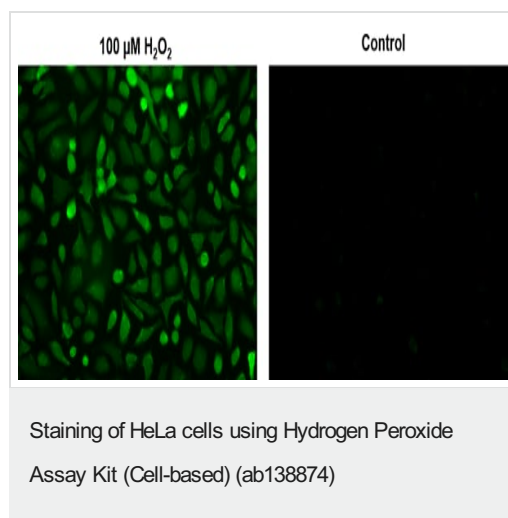
Store at -20°C. Please refer to protocols.

内容	200 tests
AbGreen Indicator	1 vial
Assay Buffer	1 x 20ml
DMSO	1 x 200µl
Hydrogen Peroxide Solution	1 x 200µl

関連性

Hydrogen Peroxide (H₂O₂) 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.

画像



Fluorescence images of intracellular hydrogen peroxide in HeLa cells using Hydrogen Peroxide Assay Kit (Cell-based) (ab138874). HeLa cells were treated with (left) or without (right) 100 µM hydrogen peroxide at 37°C for 90 minutes.

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