

Lipid Peroxidation Assay Kit (Cell-based) ab243377

★★★★★ [1 Abreviews](#) [13 References](#) [画像数 2](#)

製品の概要

製品名	Lipid Peroxidation Assay Kit (Cell-based)
検出方法	Fluorescent
製品の概要	Lipid Peroxidation Assay Kit (Cell-based) (ab243377) uses a sensitive ratiometric Lipid Peroxidation Sensor that changes its fluorescence from red to green upon peroxidation by ROS in cells, this peroxidation-dependent shift enables the ratiometric measurement of lipid peroxidation. Our kit includes H ₂ O ₂ as a positive control treatment to induce lipid peroxidation.
アプリケーション	適用あり: Flow Cyt, FM
試験プラットフォーム	Flow cytometer, Fluorescence microscope

製品の特性

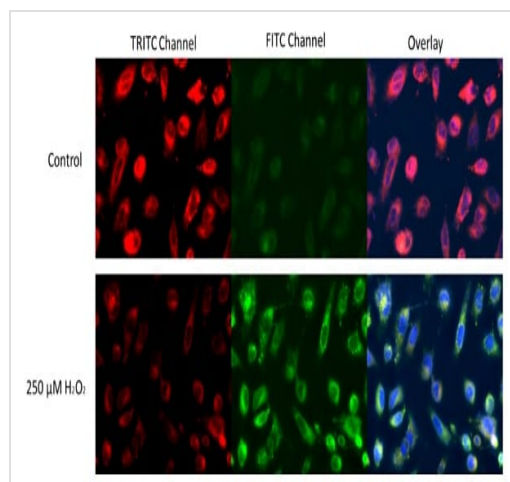
保存方法 Store at -20°C. Please refer to protocols.

内容	200 tests
3% H ₂ O ₂	1 x 100µl
HHBS	1 x 50ml
Ratiometric lipid peroxidation sensor R590/G525 (500X)	1 x 50µl

アプリケーション

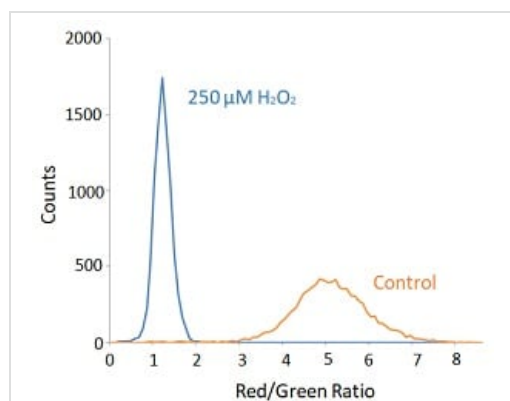
The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab243377の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Flow Cyt		Use at an assay dependent concentration.
FM		Use at an assay dependent concentration.



HeLa cells were stained with 1X Lipid Peroxidation Sensor for 30 minutes in complete growth medium at 37°C.

For H_2O_2 treatment, approximately 250 μM H_2O_2 was added to the cells and incubated for 30 minutes. The cells were then incubated with 1X Lipid peroxidation Sensor, and stained with Hoechst 33342 during the last 10 minutes of incubation. The cells were washed 3 times with HHBS and imaged with a Keyence fluorescent microscope. With H_2O_2 treatment, a clear shift of fluorescence signal of red to green was observed.



Jurkat cells were stained with 1X Lipid Peroxidation Reagent for 30 minutes in complete growth medium at 37°C.

For H_2O_2 treatment, approximately 250 μM H_2O_2 was added to the cells and incubated for 30 minutes. The cells were then incubated with 1X Lipid Peroxidation Sensor, and analyzed with a flow cytometer through FITC (488/530 nm) and PE (488/572 nm) channels. The data are represented as the ratios of red (PE)/green (FITC) fluorescence intensities. The ratio of red/green decreases in H_2O_2 treated cells indicating the presence of H_2O_2 -induced lipid peroxidation.

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