

NIR Mitochondrial Membrane Potential Assay Kit (Microplate) ab112150

★★★★★ 1 Abreviews 画像数 1

製品の概要

製品名	NIR Mitochondrial Membrane Potential Assay Kit (Microplate)
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Direct
製品の概要	<p>ab112150 is designed to detect cell apoptosis by measuring the loss of the mitochondrial membrane potential. The collapse of mitochondrial membrane potential coincides with the opening of the mitochondrial permeability transition pores, leading to the release of cytochrome C into the cytosol, which in turn triggers other downstream events in the apoptotic cascade.</p>

ab112150 NIR Mitochondria Membrane Potential Assay Kit provides all the essential components with an optimized assay method. This fluorometric assay uses our proprietary cationic MitoNIR Dye for the detection of the mitochondrial membrane potential change in cells. In normal cells, the red fluorescence intensity is increased when NIR Dye is accumulated in the mitochondria. However, in apoptotic cells, NIR stain intensity is decreased following the collapse of MMP. Cells stained with NIR Dye can be monitored fluorimetrically at $E_m = 660-680$ nm with excitation at $E_x = 620-640$ nm.

ab112150 can be used for screening apoptosis activators and inhibitors. The assay can be performed in a convenient 96-well and 384-well fluorescence microtiter-plate format.

特記事項

Related assays

Review the [cell health assay guide](#) to learn about kits to perform a [cell viability assay](#), [cytotoxicity assay](#) and [cell proliferation assay](#).

Review the [metabolism assay guide](#) to learn about assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also about how to assay metabolic function in live cells using your plate reader.

試験プラットフォーム

Microplate reader

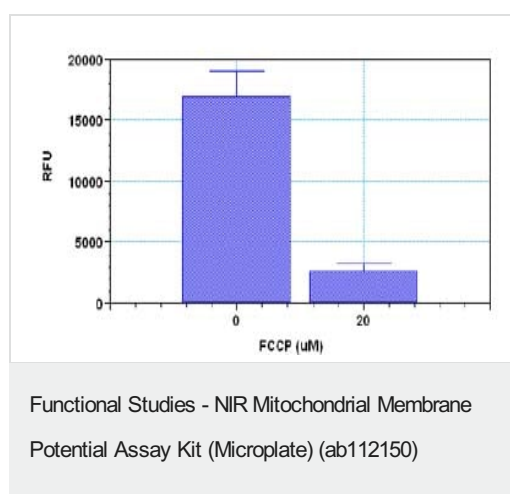
製品の特性

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

内容	5 x 96 tests
Assay Buffer A	1 x 50ml
Assay Buffer B	1 x 25ml
MitoNIR Dye	1 x 250µl

関連性 Mitochondrial Membrane Potential is an important parameter of mitochondrial function used as an indicator of cell death. The collapse of the mitochondrial Membrane potential coincides with the opening of the mitochondrial permeability transition pores, leading to the release of cytochrome c into the cytosol, which in turn triggers other downstream events in the apoptotic cascade.

画像



The decrease in NIR fluorescence with the addition of FCCP in HeLa cells.

HeLa cells were dye loaded with NIR alone, or in the presence of 20 µM FCCP for 15 minutes. The fluorescence intensity of NIR was measured 30 minutes after adding Assay buffer B (Component C) with a microplate reader at Ex/Em = 640/680 nm (cut off 665 nm).

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