

Orange Mitochondrial Membrane Potential Assay Kit (Flow Cytometry) ab138898

[1 References](#) [1 图像](#)

概述

产品名称	Orange Mitochondrial膜Potential Assay试剂盒(Flow Cytometry)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Direct
产品概述	<p>Orange Mitochondrial Membrane Potential Assay Kit (Flow Cytometry) (ab138898) is designed to detect cell apoptosis by measuring the loss of the mitochondrial membrane potential (MMP). 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>

ab138898 provides all the essential components with an optimized assay method. This fluorimetric assay uses our proprietary cationic MitoOrange Dye for the detection of apoptosis in cells with the loss of mitochondrial membrane potential. In normal cells, the red fluorescence intensity is increased when MitoOrange Dye is accumulated in the mitochondria. However, in apoptotic cells, the fluorescence intensity of MitoOrange Dye is decreased following the collapse of MMP. Cells stained with MitoOrange Dye can be visualized with a flow cytometer at 488 nm excitation with red emission (FL2 channel). The kit can be used together with other reagents for multi-parametric study of cell vitality and apoptosis. The kit is optimized for screening apoptosis activators and inhibitors with a flow cytometer.

说明

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.

平台

Fluorescence microscope

性能

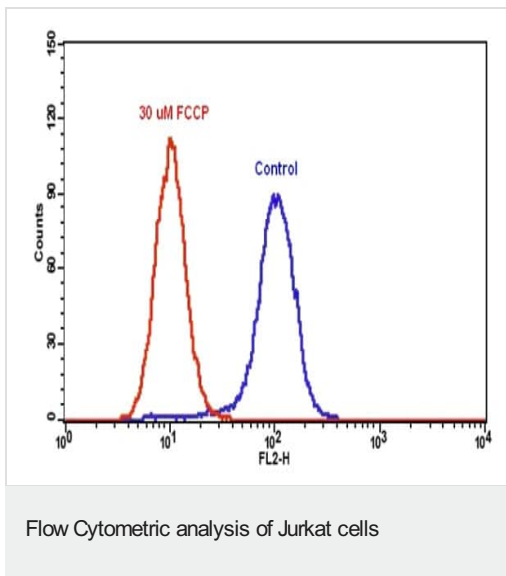
存放说明

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

组件	100 tests
Assay Buffer	1 x 100ml
MitoOrange Dye	1 x 200µl

相关性 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.

图片



Flow Cytometric analysis demonstrating the decrease in fluorescence intensity of MitoOrange Dye with the addition of FCCP in Jurkat cells. Jurkat cells were loaded with MitoOrange Dye alone (Blue) or in the presence of 30 µM FCCP (Red) for 15 minutes. The fluorescence intensity of MitoOrange Dye was measured with a flow cytometer using the FL2 channel.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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