

Mitochondrial Staining Kit - Red Fluorescence - Cytopainter ab112145

[8 References](#) [1 Image](#)

Overview

Product name	Mitochondrial Staining Kit - Red Fluorescence - Cytopainter
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Product overview	Abcam's fluorescence imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as lysosomes, mitochondria, and actin filaments. The selective staining of cell compartments provides a powerful method for studying cellular events in a spatial and temporal context.

ab112145 is designed to label mitochondria in live cells with red fluorescence. The kit uses our proprietary dye that selectively accumulates in mitochondria probably via the mitochondrial membrane potential gradient. The red fluorescent mitochondrial stain used in the kit has Ex/Em = 580/600 nm. The mitochondrial indicator, a hydrophobic compound, easily permeates intact live cells and becomes trapped in mitochondria after it gets into cells. This fluorescent mitochondrial indicator is retained in mitochondria for a long time since it carries a cell-retaining group. This key feature significantly increases the staining efficiency.

ab112145 can be readily adapted for many different types of fluorescence platforms, such as microplate assays, immunocytochemistry and flow cytometry. It is useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. The kit provides all the essential components and can be used for both proliferating and non-proliferating cells.

Review other dyes and kits for [mitochondrial staining](#), or the [live cell staining fluorescent dyes guide](#)

Notes	Mitochondria are membrane-enclosed organelles found in most eukaryotic cells. Mitochondria are sometimes described as “cellular power plants” because they generate most of the cellular supply of ATP. In addition to supplying cellular energy, mitochondria are involved in a range of other processes, such as signaling, cellular differentiation, cell death, as well as the control of the cell cycle and cell growth. Mitochondria have been implicated in several human diseases, including mitochondrial disorders and cardiac dysfunction, and may play a role in the aging process. Although most of a cellular DNA is contained in the cell nucleus, the mitochondrion has its own independent genome.
--------------	---

Platform Fluorescence microscope

Properties

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

Components	500 tests
Live Cell Staining Buffer	1 x 50ml
MitoRed Indicator	1 x 100µl

Images

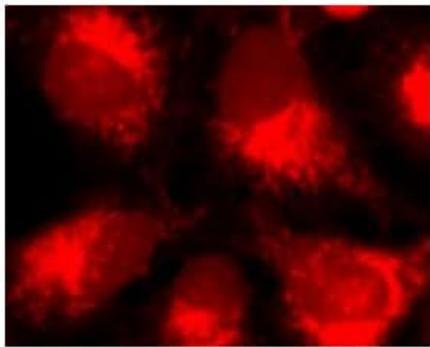


Image of U2OS cells stained with ab112145 in a black 96-well plate.

Fluorescence Microscopy- CytoPainter
Mitochondrial Staining Kit - Red Fluorescence
(ab112145)

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.com/abpromise> or contact our technical team.

Terms and conditions

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