

Mitochondrial Staining Reagent - Orange - Cytopainter ab176831

[2 Images](#)

Overview

Product name	Mitochondrial Staining Reagent - Orange - Cytopainter
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (qualitative)
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Abcam's Mitochondrial Staining Reagent - Orange Cytopainter (ab176831) is part of a series of new fluorogenic probes to label mitochondria of live cells. The proprietary mitochondria dye selectively accumulates in mitochondria probably via the mitochondrial membrane potential gradient. The mitochondrial indicator is a hydrophobic compound that easily permeates intact live cells and becomes trapped in mitochondria after it gets into cells. The 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; make it useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity.</p>

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

Notes	<p>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.</p>
--------------	--

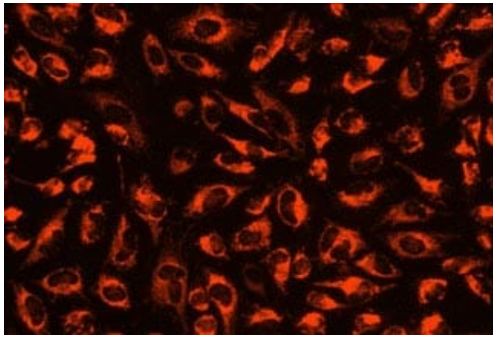
Platform	Fluorescence microscope
-----------------	-------------------------

Properties

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

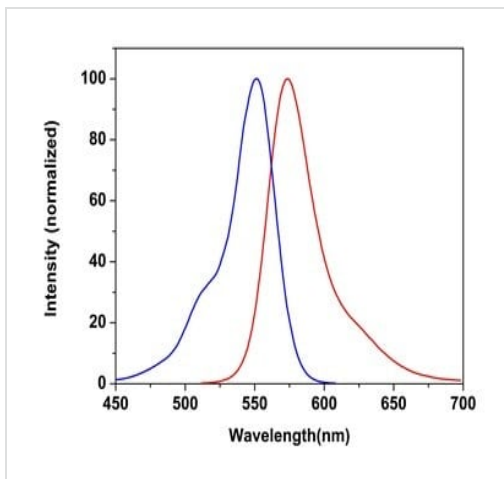
Components	500 tests
Cytopainter MitoOrange Indicator Reagent	1 x 500 tests

Images



U2OS cells stained with ab176831

Image of U2OS cells stained with Abcam's CytoPainter MitoOrange Indicator Reagent (ab176831) in a Costar black wall/clear bottom 96-well plate.



Excitation and Emission spectra for Cytopainter MitoOrange Indicator Reagent (ab176831)

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