

# Mitochondrial Staining Kit - Green Fluorescence - Cytopainter ab112143

[4 References](#) [2 Images](#)

## Overview

---

<b>Product name</b>	Mitochondrial Staining Kit - Green Fluorescence - Cytopainter
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based
<b>Product overview</b>	<p>Mitochondrial Staining Kit - Green Fluorescence   Cytopainter (ab112143) is designed to label mitochondria in live cells with green fluorescence at Ex/Em = 490/520 nm. The kit uses a proprietary dye that selectively accumulates in mitochondria probably via the mitochondrial membrane potential gradient. The mitochondrial indicator, a hydrophobic compound, easily permeates intact live cells and becomes trapped in mitochondria after it gets into cells. The 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.</p> <p>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.</p> <p>The dye will stain live cells but is not well-retained after fixation.</p> <p>Review other dyes and kits for <a href="#">mitochondrial staining</a>, or the <a href="#">live cell staining fluorescent dyes guide</a></p>
<b>Notes</b>	<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 signalling, 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>
<b>Platform</b>	Fluorescence microscope

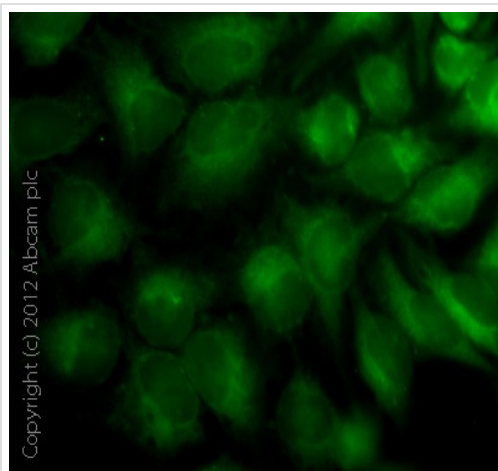
## Properties

## Storage instructions

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

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

## Images

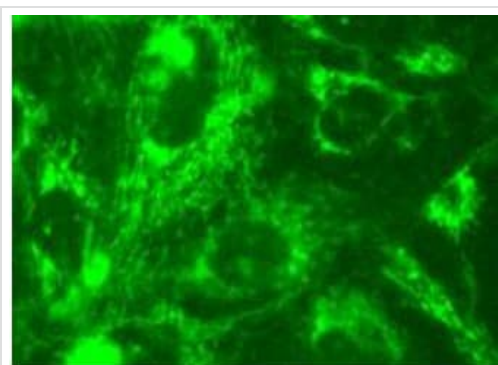


Immunocytochemistry/ Immunofluorescence -  
CytoPainter Mitochondrial Staining Kit - Green  
Fluorescence (ab112143)

### CytoPainter Mitochondrial Staining Kit - Green Fluorescence

(ab112143). HeLa cells were incubated for 60 minutes at 37C with ab112143. The staining solution was replaced with 4%PFA for 10 minutes before a final wash in PBS prior to mounting slides.

*Images obtained with an Olympus BX61 microscope using a FITC filter (495/519 nm) - 1750ms exposure.*



Immunocytochemistry/ Immunofluorescence -  
CytoPainter Mitochondrial Staining Kit - Green  
Fluorescence (ab112143)

### CytoPainter Mitochondrial Staining Kit - Green Fluorescence

(ab112143). Image of U2OS cells stained with ab112143

CytoPainter Mitochondrial Staining Kit - Green Fluorescence in a black 96-well plate.

**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