

Product datasheet

Mitochondrial Staining Kit - Blue - Cytopainter ab219939

[1 Image](#)

Overview

Product name	Mitochondrial Staining Kit - Blue - Cytopainter
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Species reactivity	Reacts with: Mammals, Other species
Product overview	Mitochondrial Staining Kit - Blue Cytopainter (ab219939) uses a proprietary fluorescent blue dye that selectively accumulates in mitochondria in live cells and can be detected at Ex/Em = 360/445 nm (DAPI filter-compatible). The mitochondrial indicator is a hydrophobic compound that easily permeates intact live cells and becomes trapped in mitochondria, where it can be retained in mitochondria for a long time thanks to its cell-retaining group. The kit can be easily adapted for many different types of fluorescence platforms, such as microplate assays, and flow cytometry.

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
-----------------	-------------------------

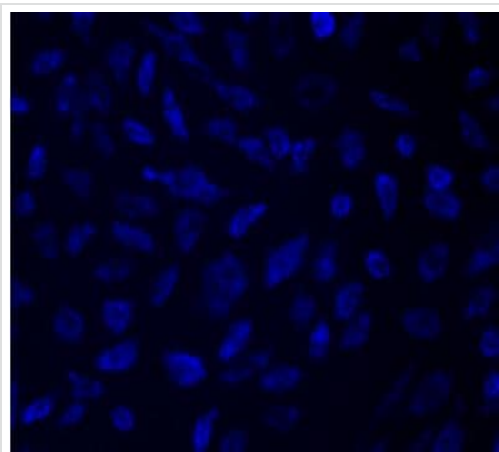
Properties

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

Components	500 tests
Live Cell Staining Buffer	1 x 50ml

Components	500 tests
MitoBlue Indicator	1 x 100µl

Images



CytoPainter Mitochondrial Staining Kit - Blue
Fluorescence (ab219939)

HeLa cells stained with CytoPainter Mitochondrial Staining Kit - Blue Fluorescence (ab219939) in a 96 well clear-bottom 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