

Product datasheet

Live Cell Labeling Kit - Blue Fluorescence - Cytopainter ab187963

[1 Image](#)

Overview

Product name	Live Cell Labeling Kit - Blue Fluorescence - Cytopainter
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Assay time	2h 00m
Product overview	<p>Live Cell Labeling Kit - Blue Fluorescence Cytopainter (ab187963) is designed to uniformly label live cells in blue fluorescence with a proprietary dye whose fluorescence is strongly enhanced upon entering into live cells. The dye is a hydrophobic compound that easily permeates intact live cells. The hydrolysis of the weakly fluorescent substrate by intracellular esterases generates a strongly fluorescent hydrophilic product that is well-retained in the cell cytoplasm. Cells grown on black wall/clear bottom plates or slides can be stained and quantified in less than two hours. View cells at Ex/Em = 360/445 nm.</p> <p>This kit can be readily adapted for a wide variety of fluorescence platforms such as microplate assays, flow cytometry and fluorescence microscope. 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 suspension and adherent cells.</p>
Notes	<p>Abcam's CytoPainter Live Cell Labeling Kits are a set of tools used to label cells for fluorescence microscopic investigations and flow cytometric investigations of cellular functions. The effective labelling of cells provides a powerful method for studying cellular events in a spatial and temporal context.</p>
Platform	Flow cytometer, Fluorescence microscope

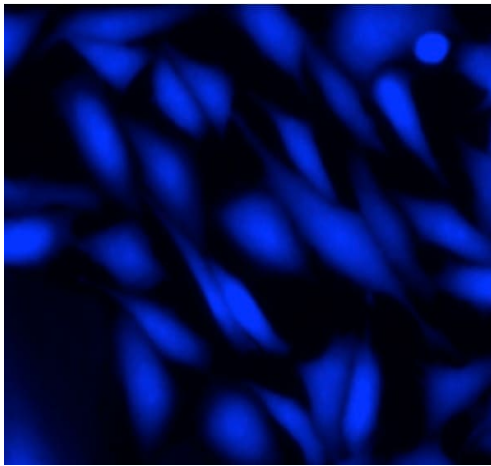
Properties

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

Components	200 tests
HHBS	1 x 100ml

Components	200 tests
Labeling Dye Blue	1 vial

Images



CPA cells

Image of CPA cells stained with CytoPainter Live Cell Labeling Kit - Blue Fluorescence (ab187963) in a black wall/clear bottom 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