

Product datasheet

Live Cell Labeling Kit - Green Fluorescence - Cytopainter ab187964

[1 Image](#)

Overview

Product name	Live Cell Labeling Kit - Green 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 - Green Fluorescence Cytopainter (ab187964) is designed to uniformly label live cells in green fluorescence. The kit uses non-fluorescent Labelling Dye Green that becomes strongly fluorescent upon entering into live cells. Labelling Dye Green is a hydrophobic compound that easily permeates intact live cells. The hydrolysis of the non-fluorescent Labelling Dye Green by intracellular esterases generates the strongly fluorescent hydrophilic calcein 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 at Ex/Em = 495/515 nm..</p> <p>This product 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 proliferating and non-proliferating, 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

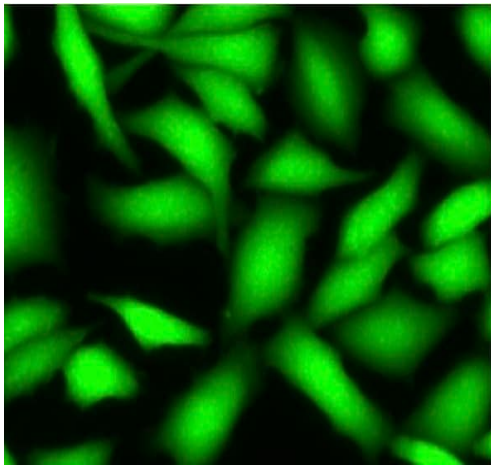
Properties

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

Components	200 tests
HHBS	1 x 100ml

Components	200 tests
Labeling Dye Green	2 vials

Images



CPA cells

Image of CPA cells stained with CytoPainter Live Cell Labeling Kit - Green Fluorescence (ab187964) 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