

Product datasheet

Live Cell Labeling Kit - Red Fluorescence - Cytopainter ab187965

[1 Image](#)

Overview

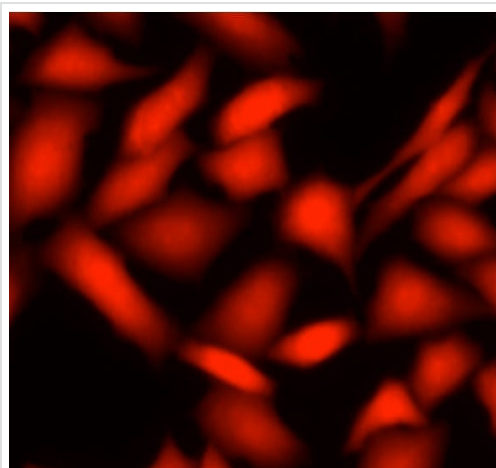
Product name	Live Cell Labeling Kit - Red Fluorescence - Cytopainter
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Assay time	2h 00m
Product overview	<p>Live Cell Labeling Kit - Red Fluorescence Cytopainter (ab187965) is designed to uniformly label live cells in red fluorescence with a proprietary non-fluorescent dye that becomes strongly fluorescence upon entering into live cells. The dye is a hydrophobic compound that easily permeates intact live cells. The hydrolysis of the non-fluorescent substrate by intracellular esterases generates a strongly red fluorescent hydrophilic product that is well-retained in the cell cytoplasm. Cells grown on slides or black wall/clear bottom plates can be stained and quantified in less than two hours.</p> <p>The assay is more robust than the tetrazolium salt one. It can be readily adapted for many different types of fluorescence platforms such as microplate assays, flow cytometry and fluorescence microscope with Ex/Em = 649/659 nm. It is also useful in a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. CytoPainter Live Cell Labeling Kit – Red Fluorescence (ab187965) provides all the essential components and can be used for both proliferating and non-proliferating 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
Labeling Dye Red	1 x 2 vials

Images



CPA cells

Image of CPA cells stained with CytoPainter Live Cell Labeling Kit - Red Fluorescence (ab187965) 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