

## Product datasheet

# Live Cell Labeling Kit - Green Fluorescence (Ex405nm) - Cytopainter ab187967

[1 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Live Cell Labeling Kit - Green Fluorescence (Ex405nm) - Cytopainter
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based
<b>Product overview</b>	<p>Live Cell Labeling Kit - Green Fluorescence (Ex405nm)   Cytopainter (ab187967) is designed to uniformly label live cells in green fluorescence for the studies that require the fluorescent tag molecules retained inside cells for a relatively longer time. The kit uses a non-fluorescent dye that carries a cell-retaining moiety. The dye becomes strongly fluorescent upon entering into live cells, and is trapped inside cells to give stable fluorescence signal that can be measured at Ex/Em = 410/500 when excited with a violet laser at Ex=405 nm. The dye is a hydrophobic compound that easily permeates intact live cells.</p> <p>The labeling process is robust, requiring minimal hands-on time. This product can be readily adapted for many different types 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 with an optimized cell-labeling protocol, and can be used for both proliferating and non-proliferating cells (either suspension or adherent cells).</p>
<b>Notes</b>	<p>Abcam's CytoPainter Cell Labeling Kits are a set of tools used to label cells for fluorescence microscopic investigations of cellular functions. The effective labeling of cells provides a powerful method for studying cellular events in a spatial and temporal context.</p>
<b>Platform</b>	Flow cytometer, Fluorescence microscope

### Properties

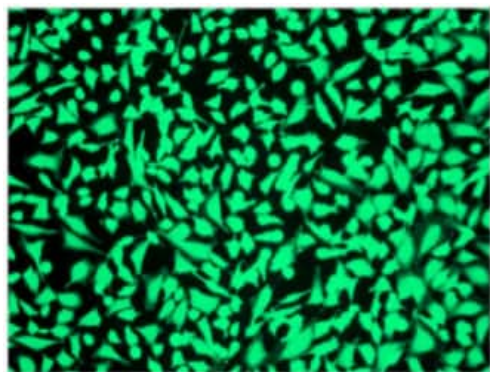
**Storage instructions** Store at -20°C. Please refer to protocols.

Components	200 tests
HHBS	1 x 100ml

Components	200 tests
Labeling Dye Green (Ex405nm)	2 vials

## Images

---



HeLa cells with CytoPainter Live Cell Labeling Kit -  
Green Fluorescence

**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