

# Cell Proliferation Staining Reagent - Green Fluorescence - Cytopainter ab176735

★★★★★ [2 Abreviews](#) [5 References](#) [画像数 2](#)

### 製品の概要

製品名	Cell Proliferation Staining Reagent - Green Fluorescence - Cytopainter
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based (quantitative)
種交差性	交差種: Mammals, Other species
製品の概要	<p>Cell Proliferation Staining Reagent - Green Fluorescence   Cytopainter (ab176735) is a powerful tool to analyze heterogeneous cell populations where multicolor analysis is required. The CytoLabeling Green Reagent is designed to label cells at Em = 511-525 nm excited with a blue laser at Ex = 488 nm.</p> <p>The CytoLabeling Green Reagent dye contained in this kit is functionally similar to CFSE and can be used for the multicolor applications where either GFP or FITC-labeled antibody is used.</p> <p>Main features of CytoLabeling Green Reagent:</p> <p><i>It eliminates the dye efflux drawback associated with CFSE</i></p> <p><i>It is compatible with the cell culture medium in the staining cells prior analysis - no washing required</i></p> <p><i>It is retained up to 9 generations after staining</i></p> <p><i>Fluorescence is preserved upon formaldehyde fixation</i></p> <p>This product can be used for imaging or flow cytometric analysis.</p> <p>Review other dyes and kits for <a href="#">cell proliferation staining</a>, or the <a href="#">live cell staining fluorescent dyes guide</a></p>
特記事項	<p>Uniform labeling of live cells with a dye containing a cell-retaining group to allow retention in proliferating cells. It enables multicolor labeling and functional analysis of live cells in combination with other fluorochromes. The dye is retained for up to 9 generations after staining.</p>

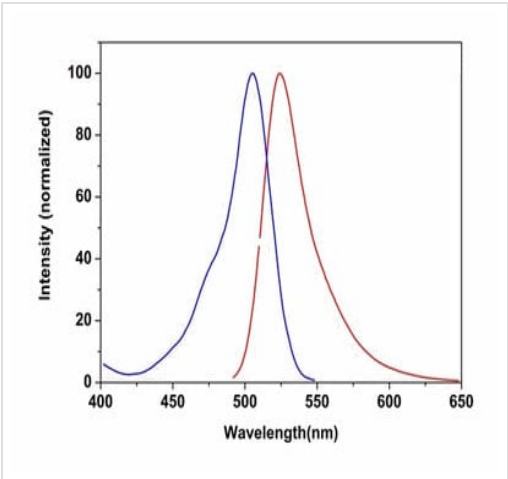
製品の特性

保存方法Store at -20°C. Please refer to protocols.

内容	500 tests	1000 tests
CytoLabeling Green Reagent	1 vial	2 vials

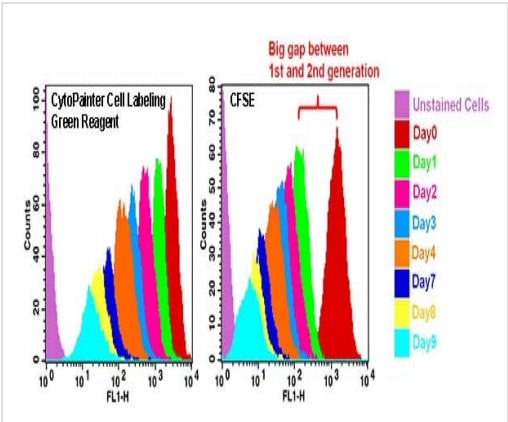
関連性Cell proliferation is the multiplication or reproduction of cells, as a result of cell growth and cell division, resulting in the expansion of a cell population.

画像



Emission and Excitation spectra for CytoPainter Cell Proliferation Staining Reagent – Green Fluorescence (ab176735).

Emission and Excitation spectra for CytoPainter Cell Proliferation Staining Reagent – Green Fluorescence (ab176735)



Cell tracking assay with CytoPainter Cell Proliferation Staining Reagent - Green Fluorescence (ab176735) and CFSE. Jurkat cells (2x10<sup>6</sup> cells/mL) were stained with ab176735 or CFSE (0.5 μM) on Day 0. The cells were passed serially at 1:1 ratio for 9 days. Fluorescence intensity was measured with a flow cytometer in FL1 channel on the day after passage. Successive generations were represented by different colors.

Cell tracking assay with CytoPainter Cell Proliferation Staining Reagent - Green Fluorescence (ab176735)

**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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors