


Product datasheet

Cell Proliferation Staining Reagent - Deep Red Fluorescence - Cytopainter ab176736

★★★★★ [2 Abreviews](#) [7 References](#) [2 Images](#)

Overview

Product name	Cell Proliferation Staining Reagent - Deep Red Fluorescence - Cytopainter
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Species reactivity	Reacts with: Other species Predicted to work with: Mammals 
Product overview	<p>Cell Proliferation Staining Reagent - Deep Red Fluorescence Cytopainter (ab176736) is a powerful tool to analyze heterogeneous cell populations where multicolor analysis is required. This reagent is designed to label cells at Em = 628-643 nm excited with a red laser at Ex = 633 nm.</p> <p>The CytoLabeling Red 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, since the CytoLabeling Reagents staining reagents have either excitation and emission spectra distinct from CFSE and its fluorescein analogs.</p> <p>Main features of CytoLabeling Deep Red 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 cell proliferation staining, or the live cell staining fluorescent dyes guide</p>
Notes	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.

Platform Flow cytometer

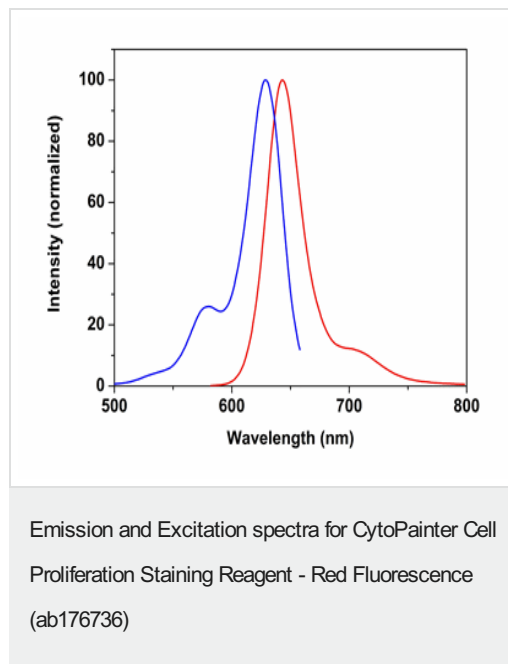
Properties

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

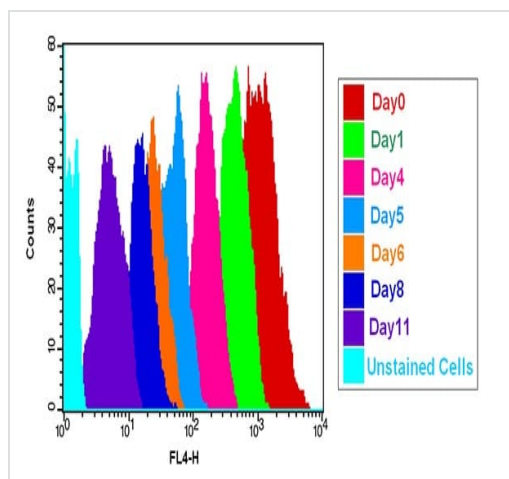
Components	500 tests	1000 tests
CytoLabeling Red Reagent	1 vial	1 x 2 vials

Relevance 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.

Images



Emission and Excitation spectra for CytoPainter Cell Proliferation Staining Reagent - Red Fluorescence (ab176736).



Cell tracking assay with CytoPainter Cell Proliferation Staining Reagent - Red Fluorescence (ab176736)

Cell tracking assay with CytoPainter Cell Proliferation Staining Reagent - Red Fluorescence (ab176736). Jurkat cells (2×10^6 cells/mL) were stained with ab176736 on Day 0. The cells were passed serially at 1:1 ratio for 11 days. Fluorescence intensity was measured with a flow cytometer in FL1 channel on the day after passage. Successive generations were represented by different colors.

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