

Product datasheet

Cell Proliferation Staining Reagent - Orange Fluorescence - Cytopainter ab176737

1 References

Overview

Product name	Cell Proliferation Staining Reagent - Orange Fluorescence - Cytopainter
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Species reactivity	Reacts with: Mammals, Other species
Product overview	Cell Proliferation Staining Reagent - Orange Fluorescence Cytopainter (ab176737) is a powerful tool to analyze heterogeneous cell populations where multicolor analysis is required. This reagent is designed to label cells at Em = 542-556 nm excited with a blue laser at Ex = 488 nm or with a green laser at Ex = 531 nm.

The CytoLabeling Orange 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.

Main features of CytoLabeling Orange Reagent:

It eliminates the dye efflux drawback associated with CFSE

It is compatible with the cell culture medium in the staining cells prior analysis - no washing required

It is retained up to 9 generations after staining

Fluorescence is preserved upon formaldehyde fixation

This product can be used for imaging or flow cytometric analysis.

Review other dyes and kits for [cell proliferation staining](#), or the [live cell staining fluorescent dyes guide](#)

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 Orange Reagent	1 vial	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.

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