

Product datasheet

Cell Tracking Dye Kit - Orange - Cytopainter ab138892

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

Overview

| | |
|-------------------------|---|
| Product name | Cell Tracking Dye Kit - Orange - Cytopainter |
| Detection method | Fluorescent |
| Sample type | Adherent cells, Suspension cells |
| Assay type | Cell-based |
| Product overview | Cell Tracking Dye Kit - Orange Cytopainter (ab138892) is designed to uniformly label live cells for studies where multicolor analysis might be required, such as cell drug resistance, cell adhesion or chemotaxis. Dye can be visualized with an orange filter (Ex/Em = 540/560 nm). |

The labelling process is robust, requiring minimal hands-on time. This product can be used in different types of fluorescence platforms such as microplate assays, flow cytometry and fluorescence microscope. The kit provides all the essential components with an optimized cell-labelling protocol, and can be used for both proliferating and non-proliferating cells (either suspension or adherent cells).

Notes Uniform labeling of live cells with a hydrophobic dye whose fluorescence is strongly enhanced upon entering the cell. Cellular esterases convert the substrate into a strong fluorescent photostable hydrophilic dye that is well retained in the cytoplasm. Cells need to be viewed in less than 24 - 48 hours as dye will be passed to daughter cells.

If you would like to track cells during several generations, we recommend **CytoPainter Cell Proliferation Staining Reagent - Orange Fluorescence (ab176737)**.

Platform Microplate reader, Fluor. microscope, Flow cyt.

Properties

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

| Components | 200 tests |
|---------------------|-----------|
| Assay Buffer | 1 x 20ml |
| Tracking dye Orange | 1 x 50µl |

Images



Fluorescence Microscopy

Image of HeLa cells stained with 1X CytoPainter Cell Tracking Staining Kit - Orange Fluorescence 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