

Product datasheet

Cell Plasma Membrane Staining Kit - Orange Fluorescence - Cytopainter ab219941

1 Image

Overview

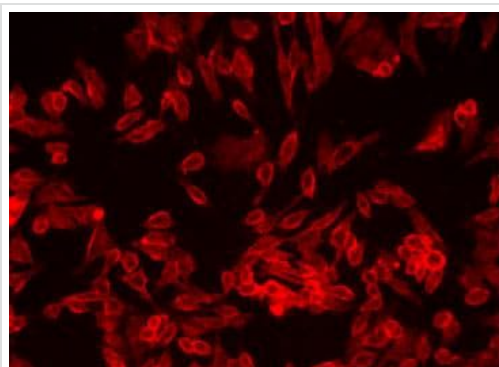
Product name	Cell Plasma Membrane Staining Kit - Orange Fluorescence - Cytopainter
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Cell Plasma Membrane Staining Kit - Orange Fluorescence Cytopainter (ab219941) uses an orange cell membrane probe (Ex/Em = 540/590 nm) that enables the uniform staining of cell membrane across a wide variety of mammalian cell types. The kit provides an excellent tool for the rapid staining of plasma membranes in suspended or attached live cells depending on the cell type and experimental conditions. The fluorescence staining in cell membranes is also maintained after fixation with formaldehyde, enabling further multi-color staining. In addition, the kit provides robust and flexible staining in live and fixed cells, and can be adapted for many different types of fluorescence platforms, such as fluorescence imaging and flow cytometry.</p> <p>Review other dyes and kits for plasma membrane staining, or the live cell staining fluorescent dyes guide</p>
Notes	<p>The cell membrane (plasma membrane) is a thin semi-permeable membrane that separates the interior of all cells from the environment. The basic function of the cell membrane is to protect the cell from its surroundings. It is composed mainly of lipids and proteins. Cell membranes are involved in a variety of cellular processes such as cell adhesion, ion conductivity and cell signaling and serve as the attachment surface for several extracellular structures, including the cell wall, glycocalyx, and intracellular cytoskeleton.</p>
Platform	Fluorescence microscope

Properties

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

Components	500 tests
Assay Buffer	1 x 50ml
DMSO	1 x 200µl
Orange Dye	1 vial

Images



CytoPainter Cell Plasma Membrane Staining Kit -
Orange Fluorescence (ab219941)

HeLa cells stained with CytoPainter Cell Plasma Membrane Staining Kit – Orange Fluorescence (ab219941) in a 96 well clear-bottom 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