

Lysosome/ER/Nuclear Staining Reagent - Cytopainter

ab139488

1 Image

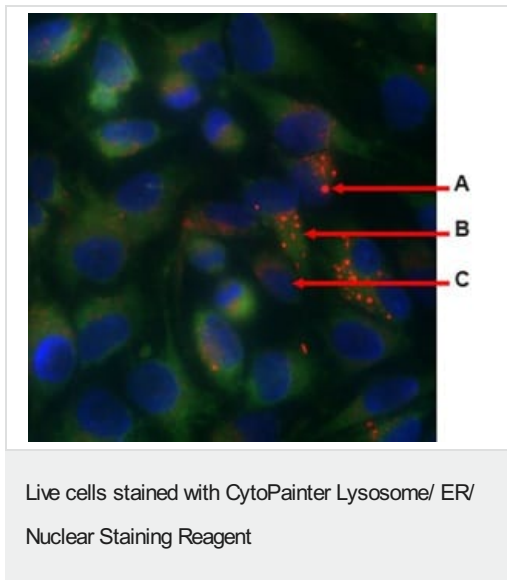
Overview

Product name	Lysosome/ER/Nuclear Staining Reagent - Cytopainter
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Assay time	0h 30m
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Lysosome/ER/Nuclear Staining Reagent Cytopainter (ab139488) containsOrganelle Reagent II, a mixture of cell-permeable red fluorescent lysosomal dye, green fluorescent endoplasmic reticulum dye and blue fluorescent nucleic acid dye. The staining pattern arising from the combination of these three dyes permits visualization of the target organelles by fluorescence/confocal microscopy. The reagent, supplied as a 500X solution, is sufficient for 1000 microscopy assays. The single-tube format makes this multi-organelle stain reagent easy to use.</p> <p>Review other dyes and kits for lysosomal staining, or the live cell staining fluorescent dyes guide</p>
Notes	<p>Lysosomal Staining - Red: Ex/Em = 568/667 nm</p> <p>Endoplasmic Reticulum Staining - Green: Ex/Em = 440/565 nm</p> <p>Nuclear Staining - Blue: Ex/Em = 350/461 nm</p>
Platform	Fluorescence microscope

Properties

Storage instructions Please refer to protocols.

Components	1000 tests
Organelle Reagent II	1 x 200µl



Live cells treated with Chloroquine and stained with Organelle Reagent II showing composite of (A) lysosomes in red, (B) endoplasmic reticula in green, and (C) nuclei in blue.

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