

Product datasheet

Lysosomal Staining Kit - Deep Red - Cytopainter
ab138896

1 References 1 Image

Overview

Product name	Lysosomal Staining Kit - Deep Red - Cytopainter
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Abcam's fluorescence imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as membranes, lysosomes, mitochondria, nuclei, etc. The selective labeling of live cell compartments provides a powerful method for studying cellular events in a spatial and temporal context.</p> <p>ab138896 is designed to label lysosomes of live cells in red fluorescence at Ex/Em = 590/620 nm. The proprietary lysotropic dye used in the kit selectively accumulates in lysosomes probably via the lysosome pH gradient. The lysotropic indicator is a hydrophobic compound that easily permeates intact live cells, and trapped in lysosomes after it gets into cells. Its fluorescence is significantly enhanced upon entering lysosomes.</p> <p>The LysoDeep Red Indicator dye used in the kit has extremely high photostability as well as excellent cellular retention makes it useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. It is suitable for proliferating and non-proliferating cells, and can be used for both suspension and adherent cells.</p> <p>Review other dyes and kits for lysosomal staining, or the live cell staining fluorescent dyes guide</p>
Notes	<p>Lysosomes are cellular organelles which contain acid hydrolase enzymes to break up waste materials and cellular debris. Lysosomes digest excess or worn-out organelles, food particles, and engulfed viruses or bacteria. The membrane around a lysosome allows the digestive enzymes to work at pH 4.5. The interior of the lysosomes is acidic (pH 4.5-4.8) compared to the slightly alkaline cytosol (pH 7.2). The lysosome maintains this pH differential by pumping protons from the cytosol across the membrane via proton pumps and chloride ion channels.</p>
Platform	Fluorescence microscope

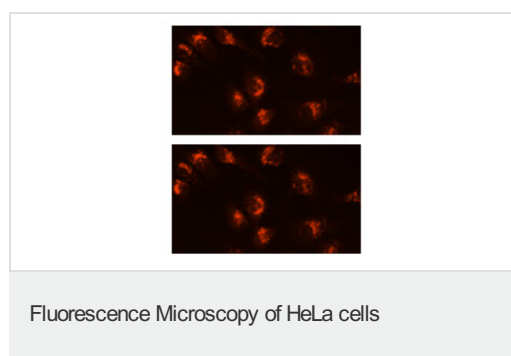
Properties

Storage instructions

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

Components	500 tests
LysoDeep Red Indicator	1 x 50µl
Staining Buffer A	1 x 25ml
Staining Buffer B	1 x 25ml

Images



HeLa cells stained with ab138896 in a black 96-well plate. The Texas Red signals were compared at 0 (upper image) and 120 seconds (lower image) exposure time by using a fluorescence microscope.

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