

Product datasheet

Lysosomal Staining Kit - Green Fluorescence (Ex405nm) - Cytopainter ab112138

1 Image

Overview

Product name	Lysosomal Staining Kit - Green Fluorescence (Ex405nm) - Cytopainter
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based
Product overview	<p>Abcam's fluorescence imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as lysosomes, mitochondria, and actin filaments. The selective staining of cell compartments provides a powerful method for studying cellular events in a spatial and temporal context.</p> <p>ab112138 is designed to label lysosomes in live cells with green fluorescence of large Stokes Shift at Ex/Em = 405/520 nm. ab112138 uses a proprietary lysotropic dye that selectively accumulates in lysosomes probably via the lysosome pH gradient. The lysotropic indicator, a hydrophobic compound, easily permeates intact live cells and gets trapped inside the lysosomes. Its fluorescence is significantly enhanced upon entering lysosomes. This key feature significantly reduces its staining background and makes it useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. The kit provides all the essential components. It is suitable for both suspension and adherent cells.</p>

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

Notes ab112138 should be stored desiccated.

Platform Fluorescence microscope

Properties

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

Components	500 tests
Live Cell Staining Buffer	1 x 50ml
LysoGreen StrK Indicator	1 x 100µl

Images

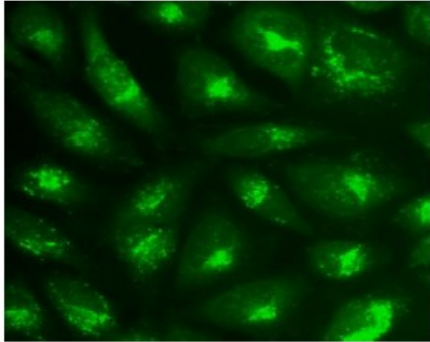


Image of U2OS cells stained with ab112138, 405 nm Excitation in a black 96-well plate

Immunocytochemistry/ Immunofluorescence -
CytoPainter Lysosomal Staining Kit - Green
Fluorescence (Ex405nm) (ab112138)

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