

Lysosomal Staining Kit - Green Fluorescence - Cytopainter

ab112136

3 References 1 图像

概述

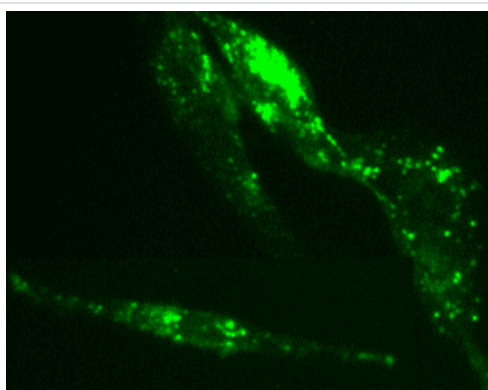
产品名称	Lysosomal Staining试剂盒- Green Fluorescence - Cytopainter
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based
产品概述	<p>Abcam's fluorescence imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as lysosomes, mitochondria, and actin filaments. The selective staining of cell compartments provides a powerful method for studying cellular events in a spatial and temporal context.</p> <p>ab112136 is designed to label lysosomes in live cells with green fluorescence at Ex/Em = 490/525 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 is trapped in lysosomes after it gets into cells. 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. 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>
平台	Fluorescence microscope

性能

存放说明 Store at -20°C. Please refer to protocols.

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

图片



ab112136 staining U2OS cells by fluorescence microscopy (in a black 96 well plate).

Fluorescence Microscopy - Lysosomal Staining Kit -
Green Fluorescence (ab112136)

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors