

Cell Navigator™ Lysosome Staining Kit *Green Fluorescence*

Catalog number: 22656 Unit size: 500 Tests

Component	Storage	Amount
Component A: LysoBrite™ Green	1 ' " " " "	1 vial (100 μL-500X DMSO stock solution)
Component B: Live Cell Staining Buffer	Freeze (<-15 °C), Minimize light exposure	1 bottle (50 mL)

OVERVIEW

Our Cell Navigator™ fluorescence imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as membranes, lysosomes, mitochondria and nuclei etc. The selective labeling of live cell compartments provides a powerful method for studying cellular events in a spatial and temporal context. This particular kit is designed to label lysosomes of live cells in green fluorescence. The kit uses a proprietary lysotropic dye that selectively accumulates in lysosomes probably vial 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. This key feature significantly increases its selectivity for lysosomes. The labeling protocol is robust, requiring minimal hands-on time. It can be readily adapted for a wide variety of fluorescence platforms such as microplate assays, immunocytochemistry and flow cytometry. It is useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. The kit provides all the essential components with an optimized cell-labeling protocol. It is suitable for proliferating and nonproliferating cells, and can be used for both suspension and adherent cells.

AT A GLANCE

Protocol summary

- 1. Prepare cells
- 2. Add LysoBrite™ Green working solution
- 3. Incubate at 37°C for 30 minutes to 2 hours
- Analyze the cells under fluorescence microscope at Ex/Em = 490/525 nm (FITC filter set)

Important Thaw all the kit components at room temperature before starting the experiment.

KEY PARAMETERS

Instrument: Fluorescence microscope

Excitation: FITC filter
Emission: FITC filter

Recommended plate: Black wall/clear bottom

PREPARATION OF WORKING SOLUTION

Add 20 µL of 500X LysoBrite™ Green stock solution (Component A) to 10 mL of Live Cell Staining Buffer (Component B) to make LysoBrite™ Green working solution. Protect from light.

Note 20 µL of 500X LysoBrite™ Green (Component A) is enough for one 96-well plate. The optimal concentration of the fluorescent lysosome indicator varies depending on the specific application. The staining conditions may be modified according to the particular cell type and the permeability of the cells or tissues to the probe.

PREPARATION OF CELL SAMPLES

For guidelines on cell sample preparation, please visit https://www.aatbio.com/resources/guides/cell-sample-preparation.html

SAMPLE EXPERIMENTAL PROTOCOL

For adherent cells:

- Grow cells either in a 96-well black wall/clear bottom plate (100 μL/well/96-well plate) or on cover-slips inside a petri dish filled with the appropriate culture medium. When cells reach the desired confluence, add equal volume of LysoBrite™ Green working solution.
- 2. Incubate the cells in a 37°C, 5% CO₂ incubator for 30 minutes to 2 hours.
- Observe the cells using a fluorescence microscope with FITC filter set (Ex/Em = 490/525 nm).

Note It is recommended to increase either the labeling concentration or the incubation time to allow the dye to accumulate if the cells do not appear to be sufficiently stained.

For suspension cells:

- Centrifuge the cells at 1000 rpm for 5 minutes to obtain a cell pellet and aspirate the supernatant.
- Resuspend the cell pellet gently in pre-warmed growth medium, and then add equal volume of LysoBrite™ Green working solution.
- 3. Incubate the cells in a 37°C, 5% $\rm CO_2$ incubator for 30 minutes to 2 hours.
- Observe the cells using a fluorescence microscope with FITC filter set (Ex/Em = 490/525 nm).

Note It is recommended to increase either the labeling concentration or the incubation time to allow the dye to accumulate if the cells do not appear to be sufficiently stained. Suspension cells may be attached to cover-slips that have been treated with BD Cell-Tak* (BD Biosciences) and stained as adherent cells.

EXAMPLE DATA ANALYSIS AND FIGURES

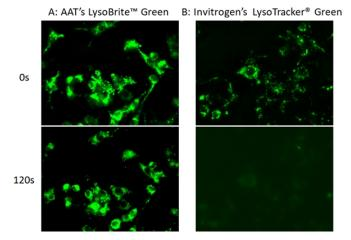


Figure 1. Images of HeLa cells stained with A: AAT's LysoBrite™ Green, B: Invitrogen's LysoTracker® Green DND-26 in a Costar black wall/clear bottom 96-well plate. Samples were continuously illuminated for 120 seconds, and the signals were compared before and after the exposure by using a Keyence fluorescence microscope.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.