

Cell Navigator™ F-Actin Labeling Kit

Orange Fluorescence

Catalog number: 22663
Unit size: 500 Tests

Component	Storage	Amount
Component A: iFluor™ 546-Phalloidin	Freeze (<-15 °C), Minimize light exposure	1 vial (50 µL)
Component B: Labeling Buffer	Freeze (<-15 °C), Minimize light exposure	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 F-actins of fixed cells in red fluorescence. The kit uses a n orange fluorescent phalloidin conjugate that is selectively bound to F-actins. This red fluorescent phalloidin conjugate is a high-affinity probe for F-actins. Used at nanomolar concentrations, phalloidins are convenient probes for labeling, identifying and quantitating F-actins in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments. The labeling protocol is robust, requiring minimal hands-on time. The kit provides all the essential components with an optimized staining protocol.

AT A GLANCE

Protocol summary

1. Prepare samples (microplate wells)
2. Remove the liquid from the plate
3. Add 100 µL/well of iFluor™ 546-Phalloidin working solution
4. Incubate the cells at room temperature for 15 to 60 minutes
5. Wash the cells
6. Examine the specimen under fluorescence microscope at Ex/Em = 550/575 nm (TRITC filter set)

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

KEY PARAMETERS

Instrument:	Fluorescence microscope
Excitation:	TRITC filter
Emission:	TRITC filter
Recommended plate:	Black wall/clear bottom

PREPARATION OF WORKING SOLUTION

Add 10 µL of iFluor™ 546-Phalloidin (Component A) to 10 mL of Labeling Buffer (Component B) to make 1X iFluor™ 546-Phalloidin working solution. Protect from light.

Note Different cell types might be stained differently. The concentration of iFluor™ 546-Phalloidin working solution should be prepared accordingly.

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

1. Perform formaldehyde fixation. Incubate the cells with 3.0% – 4.0% formaldehyde in PBS at room temperature for 10 – 30 minutes.

Note Avoid any methanol containing fixatives since methanol can disrupt

actin during the fixation process. The preferred fixative is methanol-free formaldehyde.

2. Rinse the fixed cells 2 – 3 times in PBS.
3. **Optional:** Add 0.1% Triton X-100 in PBS into fixed cells for 3 to 5 minutes to increase permeability. Rinse the cells 2 – 3 times in PBS.
4. Add 100 µL/well (96-well plate) of iFluor™ 546-Phalloidin working solution into the fixed cells.
5. Incubate the cells at room temperature for 15 to 60 minutes.
6. Rinse cells gently with PBS 2 to 3 times to remove excess dye before plate sealing.
7. Image cells using a fluorescence microscope with TRITC filter set (Ex/Em = 550/575 nm).

EXAMPLE DATA ANALYSIS AND FIGURES

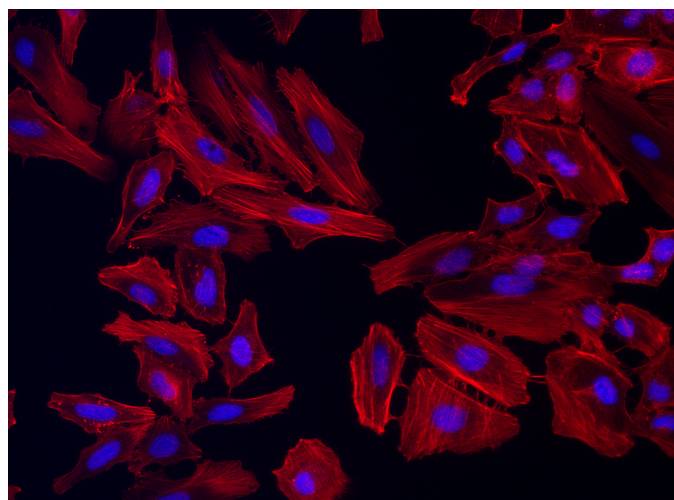


Figure 1. Fluorescence image of HeLa cells fixed with 4% formaldehyde then stained with Cell Navigator™ F-Actin Labeling Kit *Orange Fluorescence* in a Costar black 96-well plate. Cells were labeled with iFluor™ 546-Phalloidin (Cat#22663, Red) and nuclei stain DAPI (Cat#17507, Blue), respectively.

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.