

Phalloidin-iFluor® 405 Conjugate

 Catalog number: 23111
 Unit size: 300 Tests

Component	Storage	Amount (Cat No. 23111)
Phalloidin-iFluor® 405 Conjugate	Freeze (< -15 °C), Minimize light exposure	300 Tests

OVERVIEW

This blue fluorescent phalloidin conjugate (equivalent to Alexa Fluor® 405-labeled phalloidin) selectively binds to F-actins. Used at nanomolar concentrations, phalloidin derivatives are convenient probes for labeling, identifying and quantitating F-actins in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments. Phalloidin binds to actin filaments much more tightly than to actin monomers, leading to a decrease in the rate constant for the dissociation of actin subunits from filament ends, essentially stabilizing actin filaments through the prevention of filament depolymerization. Moreover, phalloidin is found to inhibit the ATP hydrolysis activity of F-actin. Phalloidin functions differently at various concentrations in cells. When introduced into the cytoplasm at low concentrations, phalloidin recruits the less polymerized forms of cytoplasmic actin as well as filamin into stable "islands" of aggregated actin polymers, yet it does not interfere with stress fibers, i.e. thick bundles of microfilaments. The property of phalloidin is a useful tool for investigating the distribution of F-actin in cells by labeling phalloidin with fluorescent analogs and using them to stain actin filaments for light microscopy. Fluorescent derivatives of phalloidin have turned out to be enormously useful in localizing actin filaments in living or fixed cells as well as for visualizing individual actin filaments in vitro. Fluorescent phalloidin derivatives have been used as an important tool in the study of actin networks at high resolution. AAT Bioquest offers a variety of fluorescent phalloidin derivatives with different colors for multicolor imaging applications.

AT A GLANCE
Protocol Summary

1. Prepare samples in microplate wells
2. Remove liquid from samples in the plate
3. Add Phalloidin-iFluor® 405 Conjugate solution (100 µL/well)
4. Stain the cells at room temperature for 20 to 90 minutes
5. Wash the cells
6. Examine the specimen under microscope with DAPI filter

Important note

Warm the vial to room temperature and centrifuge briefly before opening.

Storage and Handling Conditions

The solution should be stable for at least 6 months if store at -20 °C. Protect the fluorescent conjugates from light, and avoid freeze/thaw cycles.

Note Phalloidin is toxic, although the amount of toxin present in a vial could be lethal only to a mosquito (LD50 of phalloidin = 2 mg/kg), it should be handled with care.

PREPARATION OF WORKING SOLUTION
Phalloidin-iFluor® 405 Conjugate working solution

Add 1 µL of Phalloidin-iFluor® 405 Conjugate solution to 1 mL of PBS with 1% BSA.

Note The stock solution of phalloidin conjugate should be aliquoted and stored at -20 °C, protected from light.

Note Different cell types might be stained differently. The

concentration of phalloidin conjugate working solution should be prepared accordingly.

SAMPLE EXPERIMENTAL PROTOCOL
Stain the cells

1. Perform formaldehyde fixation. Incubate 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 Phalloidin-iFluor® 405 Conjugate working solution into the fixed cells, and stain the cells at room temperature for 20 to 90 minutes.
5. Rinse cells gently with PBS 2 to 3 times to remove excess phalloidin conjugate before plating, sealing and imaging under microscope with DAPI filter set.

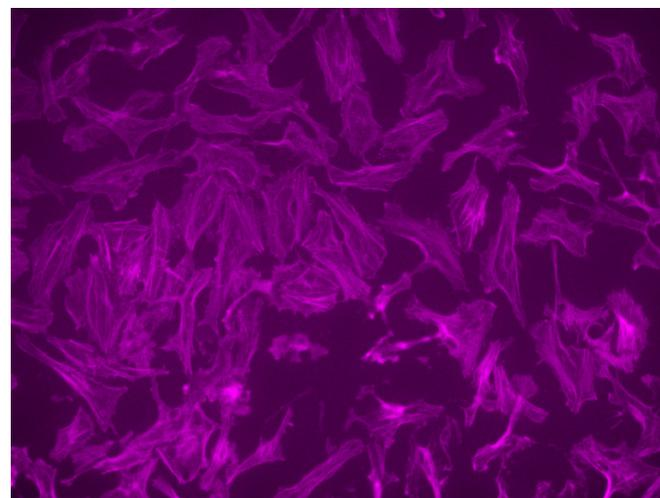
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. Fluorescence image of HeLa cells stained with Phalloidin-iFluor® 405 Conjugate using fluorescence microscope with a Violet filter set. The cells were fixed in 4% formaldehyde, permeabilised with 0.2% Triton X-100 followed by staining with Phalloidin-iFluor® 405 Conjugate. The image was acquired with fluorescence microscope using Violet filter.

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.