

**XFD647 Phalloidin**

 Catalog number: 23159  
 Unit size: 300 Tests

Component	Storage	Amount (Cat No. 23159)
XFD647 Phalloidin *equivalent to Alexa Fluor® 647 phalloidin*	Freeze (< -15 °C), Minimize light exposure	300 Tests

**OVERVIEW**

XFD647 phalloidin is a high-affinity probe for filamentous actin (F-actin), conjugated to XFD647, a bright, photostable, far-red fluorescent dye. XFD647, manufactured by AAT Bioquest, is structurally similar to Alexa Fluor™ 647 (Thermo Fisher), ensuring comparable performance in fluorescence applications. Phalloidin, a bicyclic peptide toxin derived from \*Amanita phalloides\* (commonly known as the death cap mushroom), is well known for its high specificity and strong binding affinity to F-actin. When conjugated to XFD647, phalloidin enables precise, high-contrast visualization of F-actin structures with minimal background interference. This fluorescent conjugate is highly effective for imaging F-actin in a wide range of biological specimens, including plant and animal tissues, fixed and permeabilized cells, and cell-free systems. XFD647 phalloidin binds to F-actin with nanomolar affinity, making it an excellent tool for labeling, identifying, and quantifying actin filaments. Its superior photostability and blinking properties make it particularly well suited for super-resolution microscopy techniques such as SIM and STORM. Additionally, XFD647 phalloidin staining is fully compatible with other fluorescent labels, including fluorescent proteins, Qdot nanocrystals, and iFluor®-conjugated secondary antibodies, allowing seamless integration into multiplex fluorescence imaging workflows.

**AT A GLANCE**
**Protocol Summary**

1. Prepare samples in microplate wells
2. Remove liquid from samples in the plate
3. Add XFD647 Phalloidin 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 Cy3 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.

**KEY PARAMETERS**
**Fluorescence microscope**

Emission	Cy5 filter
Excitation	Cy5 filter
Recommended plate	Black wall/clear bottom

**PREPARATION OF STOCK SOLUTIONS**

*Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles*

**XFD647 Phalloidin stock solution**

1. Add 30 µL of DMSO to the powder and mix well.
2. Add 1 µL of the XFD647 Phalloidin 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 the phalloidin conjugate working solution should be prepared accordingly.

**PREPARATION OF WORKING SOLUTION**
**XFD647 Phalloidin Conjugate working solution**

1. Add 1 µL of XFD647 Phalloidin 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 XFD647 Phalloidin 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 Cy5 filter set.

**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](mailto:info@aatbio.com) if you have any questions.