

Phalloidin-iFluor 647 Reagent ab176759

★★★★★ [5 Abreviews](#) [76 References](#) [4 图像](#)

概述

| | |
|------|--|
| 产品名称 | Phalloidin-iFluor 647试剂 |
| 样品类型 | Adherent cells, Suspension cells |
| 检测类型 | Cell-based (qualitative) |
| 产品概述 | <p>Phalloidin-iFluor 647 Reagent (ab176759) is one of a series of phalloidin conjugates that bind to actin filaments, also known as F-actin. Phalloidin-iFluor 647 can be detected with a fluorescent microscope at Ex/Em = 650/665 nm.</p> <p>Phalloidin conjugates are convenient probes for labeling, identifying and quantifying animal or plant actin filaments in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments. They can also be used in paraffin-embedded samples that have been de-paraffinized.</p> |

Review other popular phalloidin dye conjugates, including [Phalloidin-iFluor 488](#), [Phalloidin-iFluor 594](#), [Phalloidin-iFluor 555](#), and [Rhodamine Phalloidin](#), search the website to see [all phalloidin conjugates](#), or read the [phalloidin staining protocol](#).

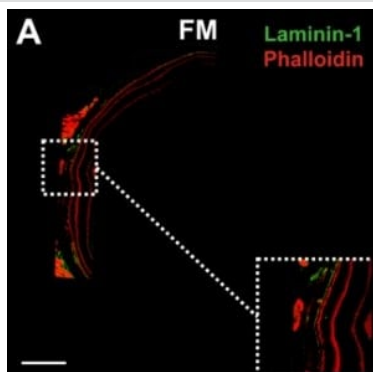
| | |
|----|---|
| 说明 | <p>Staining fixed cell or tissue samples with phalloidin conjugates is very simple; it requires a single 20-90 min incubation with the phalloidin, followed by 3 short wash steps. Phalloidin staining can be combined with antibody-based staining by adding the phalloidin conjugate during either the primary or secondary antibody incubation step.</p> <p>When used in unfixed samples, phalloidin binding leads to a decrease in the disassociation rate of actin subunits from the ends of actin filaments, essentially stabilizing actin filaments through the prevention of filament depolymerisation.</p> |
|----|---|

| | |
|----|-------------------------|
| 平台 | Fluorescence microscope |
|----|-------------------------|

性能

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

| 组件 | 300 tests |
|---------------------------------|---------------|
| Phalloidin-iFluor 647 Conjugate | 1 x 300 tests |

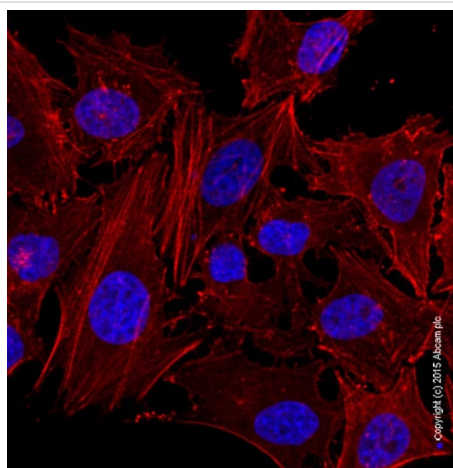


Burgoyne, Thomas et al.
PloS one vol. 13,1 e0191048.
(2018)

Functional Studies - Phalloidin-iFluor 647 Reagent
(ab176759)

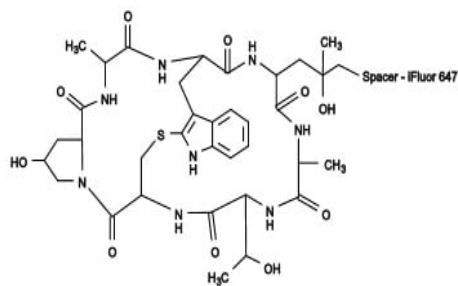
Burgoyne, Thomas et al., PloS one?vol. 13,1
e0191048., supplementary Fig 2,
doi:10.1371/journal.pone.0191048

Fluorescence microscopy (FM) image of laminin-1 and actin (phalloidin) staining.

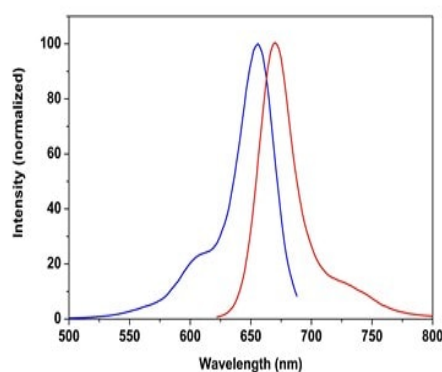


CytoPainter Phalloidin-iFluor 647 Reagent
(ab176759)

Actin filaments staining in HeLa cells. Actin filaments (red) were stained with CytoPainter Phalloidin-iFluor 647 reagent (ab176759). Nuclei were stained with DAPI (blue).



Structure



Cytopainter Phalloidin-iFluor 647 Reagent
(ab176759)

Excitation and emission spectra of phalloidin-iFluor 647 reagent.

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