

Phalloidin-iFluor 594 Reagent ab176757

★★★★★ [6 Abreviews](#) [72 References](#) [4 图像](#)

概述

产品名称	Phalloidin-iFluor 594试剂
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based (qualitative)
产品概述	<p>Phalloidin-iFluor 594 Reagent (ab176757) is one of a series of phalloidin conjugates that bind to actin filaments, also known as F-actin. Phalloidin-Fluor 594 can be detected with a fluorescent microscope at Ex/Em = 590/618 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>

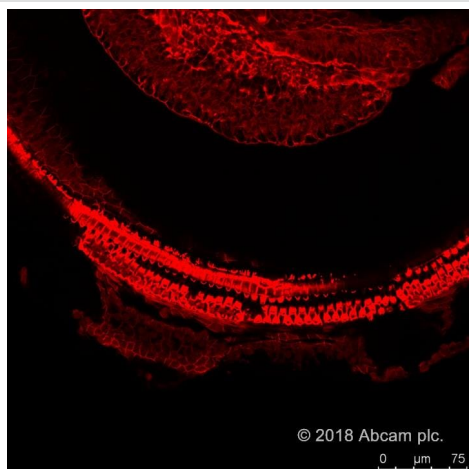
说明	<p>Review other popular phalloidin dye conjugates, including Phalloidin-iFluor 488, Phalloidin-iFluor 647, Phalloidin-iFluor 555, and Rhodamine Phalloidin, search the website to see all phalloidin conjugates, or read the phalloidin staining protocol.</p> <p>Staining fixed cell or tissue samples with 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.
------	--

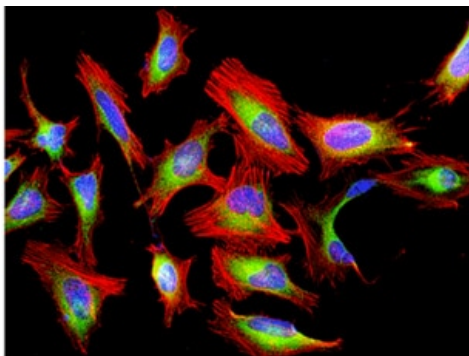
组件	300 tests
Phalloidin-iFluor 594 Conjugate	1 x 300 tests



Whole Mount of Mouse Inner Ear stained with
Phalloidin 594

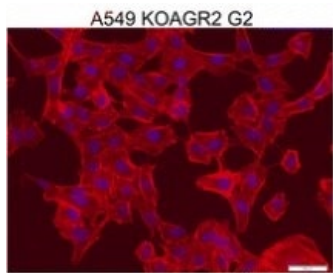
Image courtesy of Mr. Shahar Taiber

Ab176757 at 1/1000 dilution in PBS used for a whole-mount of the sensory epithelium in a mouse inner ear.



CytoPainter Phalloidin-iFluor 594 Reagent
(ab176757)

Actin filaments staining in HeLa cells. Actin filaments (red) were stained with CytoPainter Phalloidin-iFluor 594 reagent (ab176757); ER (green) was stained with CytoPainter ER Staining Kit – Green Fluorescence ([ab139481](#)). Nuclei (blue) were stained with DAPI.



F-actin/ DAPI

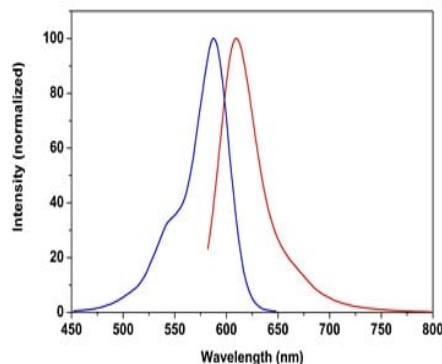
Sommerova, Lucia et al.
BMC cancer vol. 17,1
546. (2017)

The A549 KOAGR2 were stained with CytoPainter phalloidin-iFluor 594 reagent to detect morphological changes associated with F-actin reorganization. Nuclei were visualised with DAPI.

Functional Studies - Phalloidin-iFluor 594 Reagent

(ab176757)

Sommerova, Lucia et al., BMC cancer vol. 17,1 546., Fig 3, doi:10.1186/s12885-017-3537-5



Excitation and emission spectra of phalloidin-iFluor 594 reagent.

CytoPainter Phalloidin-iFluor 594 Reagent

(ab176757)

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