

Phalloidin-iFluor 555 Reagent ab176756

★★★★★ [3 Abreviews](#) [58 References](#) [3 图像](#)

概述

产品名称 Phalloidin-iFluor 555试剂

样品类型 Adherent cells, Suspension cells

检测类型 Cell-based (qualitative)

产品概述 Phalloidin-iFluor 555 Reagent (ab176756) is one of a series of phalloidin conjugates that bind to actin filaments, also known as F-actin. Phalloidin-iFluor 555 can be detected with a fluorescent microscope at Ex/Em = 556/574 nm.

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 deparaffinized.

Review other popular phalloidin dye conjugates, including [Phalloidin-iFluor 488](#), [Phalloidin-iFluor 647](#), [Phalloidin-iFluor 594](#), and [Rhodamine Phalloidin](#), search the website to see [all phalloidin conjugates](#), or read the [phalloidin staining protocol](#).

说明 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.

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.

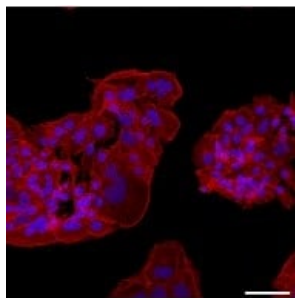
平台 Fluorescence microscope

性能

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

组件	300 tests
Phalloidin-iFluor 555 Conjugate	1 x 300 tests

图片

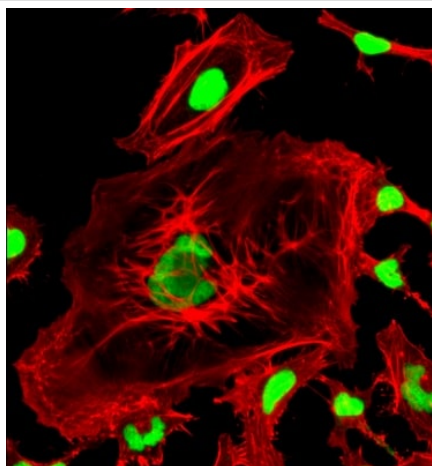


Wong, Michael K et al.
PloS one vol. 13,6
e0199632. (2018)

Functional Studies - Phalloidin-iFluor 555 Reagent
(ab176756)

Wong, Michael K et al., PloS one?vol. 13,6 e0199632.,
Fig 6, doi:10.1371/journal.pone.0199632

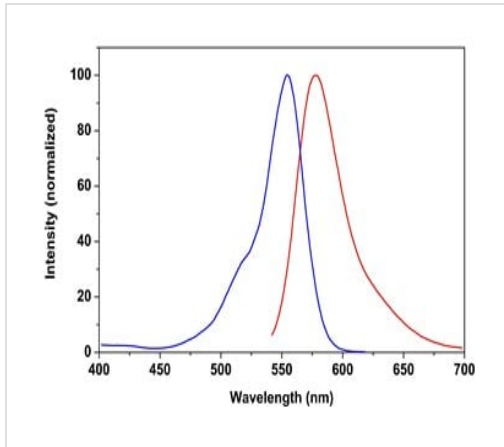
Immunofluorescent images of Phalloidin staining at day 3 across various surface thicknesses. Red fluorescence indicates phalloidin staining for F-actin and blue fluorescence indicates DAPI staining for cell nuclei.



CytoPainter Phalloidin-iFluor 555 Reagent
(ab176756)

Actin filaments staining in HeLa cells. Actin filaments (red) were stained with CytoPainter Phalloidin-iFluor 555 reagent ([ab176755](#)). Nuclei were stained with **Nuclear Green DCS1** ([ab138905](#)).

Excitation and emission of phalloidin-iFluor 555 reagent.



CytoPainter Phalloidin-iFluor 555 Reagent
(ab176756)

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