

# Phalloidin-iFluor 488 Reagent ab176753

★★★★★ [14 Abreviews](#) [215 References](#) [画像数 4](#)

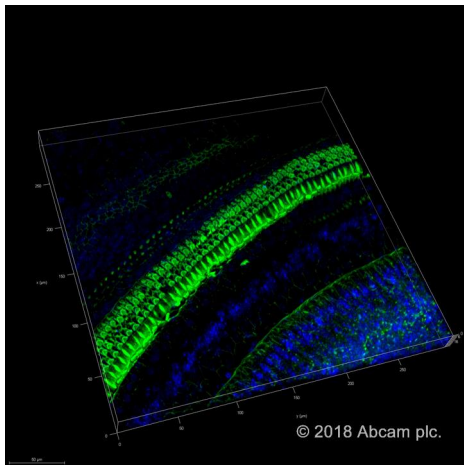
### 製品の概要

製品名	Phalloidin-iFluor 488 Reagent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based (qualitative)
製品の概要	<p>Phalloidin-iFluor 488 Reagent ab176753 is one of a series of phalloidin conjugates that bind to actin filaments, also known as F-actin. The iFluor 488 dye can be easily detected with a fluorescent microscope at Ex/Em = 493/517 nm.</p> <p>Our 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 with paraffin-embedded samples that have been de-paraffinized.</p> <p>Review other popular phalloidin dye conjugates, including <a href="#">Phalloidin-iFluor 647</a>, <a href="#">Phalloidin-iFluor 594</a>, <a href="#">Phalloidin-iFluor 555</a>, and <a href="#">Rhodamine Phalloidin</a>, search the website to see <a href="#">all phalloidin conjugates</a>, or read the <a href="#">phalloidin staining protocol</a>.</p>
特記事項	<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>
試験プラットフォーム	Fluorescence microscope

### 製品の特性

**保存方法** Store at -20°C. Please refer to protocols.

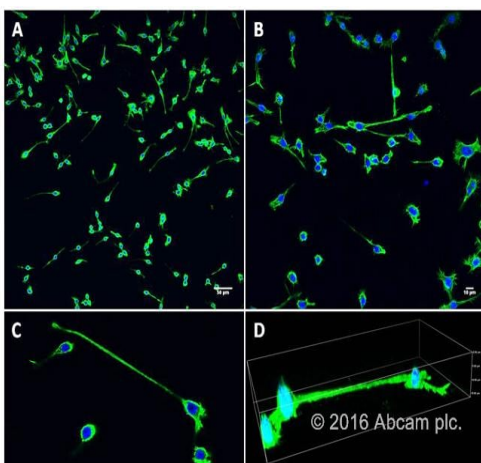
内容	300 tests
Phalloidin-iFluor 488 Conjugate	1 x 300 tests



Ab176753 at 1/1000 dilution in PBS whole mount  
Immunofluorescence of Mouse inner ear sensory epithelia. Tissue was incubated for 2 hours at room temperature.

Whole Mount Immunofluorescence of Mouse Inner Ear

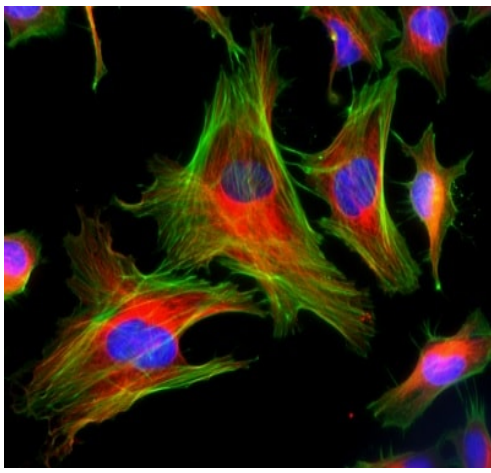
Image courtesy of Mr. Shahar Taiber



Macrophage J774A.1 cell line stained with Phalloidin 488 and DAPI nuclear counterstaining.

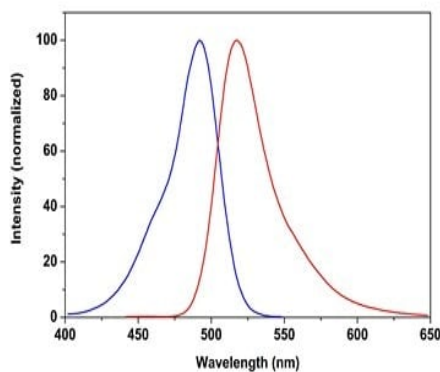
Phalloidin iFluor 488 staining of macrophages

Review by Dr. Jose Ramos Vivas



CytoPainter Phalloidin-iFluor 488 Reagent  
(ab176753)

Actin filaments staining in HeLa cells. Actin filaments (green) were stained with CytoPainter Phalloidin-iFluor 488 reagent (ab176753); tubulin filaments were stained with a mouse anti-tubulin antibody/goat anti-mouse IgG (red). Nuclei were stained with Hoechst 33342.



CytoPainter Phalloidin-iFluor 488 Reagent  
(ab176753)

Excitation and emission spectra of phalloidin-iFluor 488 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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors