

## Product datasheet

# Phalloidin-iFluor 532 Reagent ab176755

[2 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Phalloidin-iFluor 532 Reagent
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based (qualitative)
<b>Product overview</b>	<p>Phalloidin-iFluor 532 Reagent (ab176755) is one of a series of phalloidin conjugates that bind to actin filaments, also known as F-actin. Phalloidin-iFluor 532 dye can be easily detected with a fluorescent microscope at Ex/Em = 542/558 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 deparaffinized.</p> <p>Phalloidin binds to actin filaments much more tightly than to actin monomers, leading to a decrease in the dissociation rate of actin subunits from filaments ends, essentially stabilizing actin filaments through the prevention of filament depolymerization.</p>
<b>Platform</b>	Fluorescence microscope

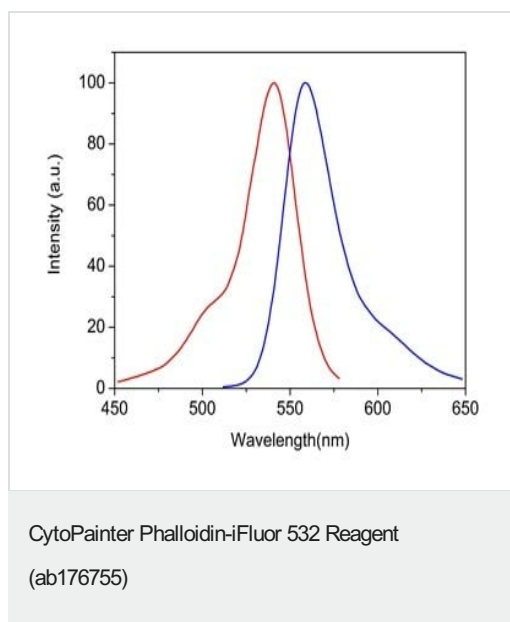
### Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	300 tests
Phalloidin-iFluor 532 Conjugate	1 x 300 tests

### Images

Excitation and emission spectra of phalloidin-iFluor 532 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.com/abpromise> or contact our technical team.

### Terms and conditions

---

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