

Phalloidin-iFluor 514 Reagent ab176754

[1 Image](#)

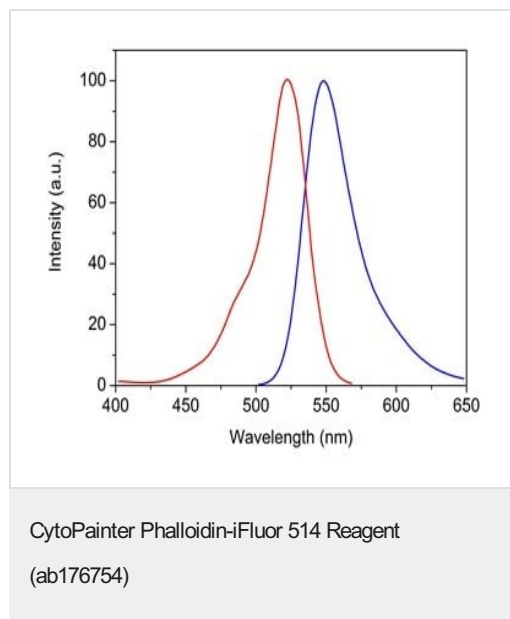
Overview

Product name	Phalloidin-iFluor 514 Reagent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (qualitative)
Product overview	<p>Phalloidin-iFluor 514 Reagent (ab176754) is one of a series of phalloidin conjugates that bind to actin filaments, also known as F-actin. Phalloidin-iFluor 514 can be easily detected with a fluorescent microscope at Ex/Em = 520/547 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> <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>
Notes	<p>Actin is a globular, roughly 42-kDa protein found in almost all eukaryotic cells. It is also one of the most highly conserved proteins, differing by no more than 20% in species as diverse as algae and humans. Actin is the monomeric subunit of two types of filaments in cells: microfilaments, one of the three major components of the cytoskeleton, and thin filaments, part of the contractile apparatus in muscle cells. Thus, actin participates in many important cellular processes including muscle contraction, cell motility, cell division and cytokinesis, vesicle and organelle movement, cell signaling, as well as the establishment and maintenance of cell junctions and cell shape.</p>
Platform	Fluorescence microscope

Properties

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

Components	300 tests
Phalloidin-iFluor 514 Conjugate	1 x 300 tests



Excitation and emission spectra of phalloidin-iFluor 514 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