

## Product datasheet

# F-actin Staining Kit - Orange Fluorescence - Cytointer ab112126

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

### Overview

<b>Product name</b>	F-actin Staining Kit - Orange Fluorescence - Cytointer
<b>Sample type</b>	Tissue, Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based
<b>Product overview</b>	<p>F-actin Staining Kit - Orange Fluorescence   Cytointer (ab112126) fluorescence imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as lysosomes, mitochondria, and actin filaments. The selective staining of cell compartments provides a powerful method for studying cellular events in a spatial and temporal context.</p> <p>ab112126 is designed to stain F-actins in fixed cells with orange fluorescence. ab112126 uses an orange fluorescent phalloidin conjugate that is selectively bound to F-actins. The orange fluorescent phalloidin conjugate has Ex/Em = 550/575 nm. The conjugate also has good photostability for fluorescence imaging. It is a high-affinity probe for labeling, identifying and quantitating F-actins in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments. ab112126 provides all the essential components with an optimized labeling protocol.</p>
<b>Notes</b>	ab112126 should be stored dessicated.
<b>Platform</b>	Fluorescence microscope

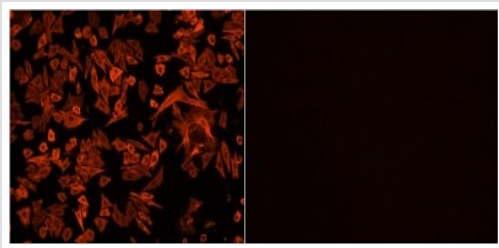
### Properties

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

Components	500 tests
Labeling Buffer	1 x 50ml
Orange Fluorescent Phalloidin Conjugate	1 x 50µl

## Images

---



Functional Studies - CytoPainter F-actin Staining Kit  
- Orange Fluorescence (ab112126)

ab112126, staining CPA cells fixed with formaldehyde in a black 96-well plate.

Left panel: The cells were labeled with Orange Fluorescent Phalloidin Conjugate for 30 minutes only.

Right panel: The cells were treated with phalloidin for 10 minutes, then stained with Orange Fluorescent Phalloidin Conjugate for 30 minutes.

**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