

F-actin Staining Kit - Orange Fluorescence - Cytopainter ab112126

1 [References](#) 1 [图像](#)

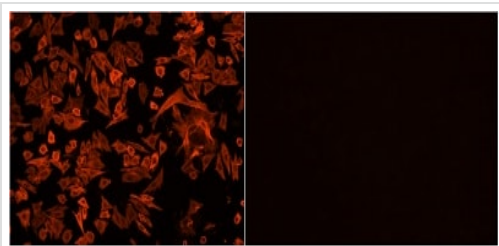
概述

产品名称	F-actin Staining试剂盒- Orange Fluorescence - Cytopainter
样品类型	Tissue, Adherent cells, Suspension cells
检测类型	Cell-based
产品概述	<p>F-actin Staining Kit - Orange Fluorescence Cytopainter (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>
说明	ab112126 should be stored dessicated.
平台	Fluorescence microscope

性能

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

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



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.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