

# F-actin Staining Kit - Green Fluorescence - Cytopainter ab112125

★★★★☆ 1 Abreviews 10 References 画像数 1

製品の概要

製品名	F-actin Staining Kit - Green Fluorescence - Cytopainter
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Cell-based (qualitative)
製品の概要	<p>F-actin Staining Kit - Green Fluorescence   Cytopainter (ab112125) 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>ab112125 is designed to stain F-actins in fixed cells with green fluorescence. The kit uses a green fluorescent phalloidin conjugate that is selectively bound to F-actins. This green fluorescent phalloidin conjugate is a high-affinity probe for F-actins. When used at nanomolar concentrations, phalloidins are convenient probes for labeling, identifying and quantitating F-actins in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments.</p> <p>ab112125 provides all the essential components with an optimized staining protocol, which is robust requiring minimal hands-on time. It is an excellent tool for preserving the fluorescent images of particular cells. The phalloidin conjugate has spectral properties similar to those of FITC (Ex/Em = 500/520 nm). It can also be used for fluorescence microscope demonstrations using FITC filter (Ex/Em = 490/525 nm).</p>
特記事項	ab112125 should be stored dessicated.
試験プラットフォーム	Fluorescence microscope

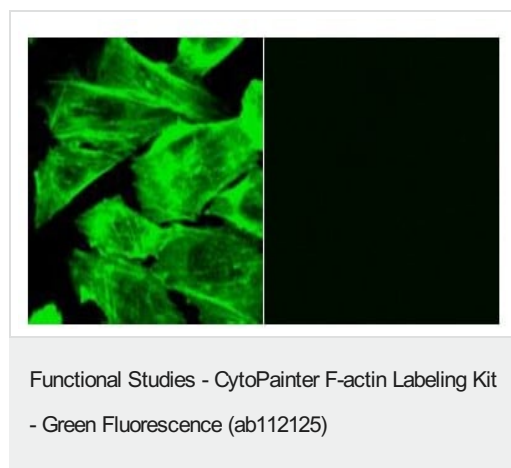
製品の特性

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

内容	500 tests

内容	500 tests
Green Fluorescent Phalloidin Conjugate	1 x 50µl
Labeling Buffer	1 x 50ml

## 画像



Left Panel: Cells labeled with 1X Green Fluorescent Phalloidin Conjugate for 30 minutes only.

Right Panel: Cells treated with phalloidin for 10 minutes, then stained with Green 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.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