

ab112127

CytoPainter F-actin Staining Kit - Red Fluorescence

Instructions for Use

For labelling sub-cellular organelles in tissue sections, cell cultures and cell free experiments using our proprietary Red fluorescence probe

This product is for research use only and is not intended for diagnostic use.

Version: 4 Last Updated: 09 December 2013

Table of Contents

1.	Introduction	3
2.	Protocol Summary	5
3.	Kit Contents	6
4.	Storage and Handling	6
5.	Assay Protocol	7
6.	Data Analysis	12

1. Introduction

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.

Abcam CytoPainter imaging kits are a set of fluorescence imaging tools for labeling sub-cellular organelles such as membranes, lysosomes, mitochondria, nuclei, etc. The selective labeling of live cell compartments provides a powerful method for studying cellular events in a spatial and temporal context.

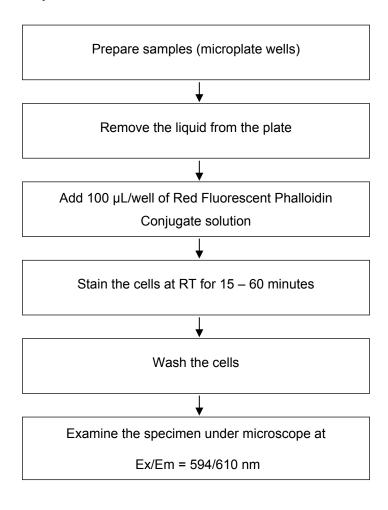
ab112127 is designed to label F-actins in fixed cells with red fluorescence. The red fluorescent phalloidin conjugate that selectively binds to F-actins. When used at nanomolar concentrations, phallotoxins are convenient probes for labeling, identifying and quantitating F-actins in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments.

Ab112127 provides all the essential components with an optimized labeling protocol.

The phalloidin conjugate has Ex/Em = 594/610 nm, compatible with Texas Red filter.

2. Protocol Summary

Summary for One 96-well Plate



Note: Warm all the components to room temperature before opening.

3. Kit Contents

Components	Amount
Component A: Red Fluorescent Phalloidin Conjugate	1 x vial
Component B: Labeling Buffer	50 mL

4. Storage and Handling

Keep at -20°C. Avoid exposure to light.

5. Assay Protocol

Note: This protocol is for one 96 - well plate.

F-ACTIN STAINING ONLY

A. Preparation of 1X Red Fluorescent Phalloidin Conjugate Working Solution

Add 10 μ L Red Fluorescent Phalloidin Conjugate (Component A) to 10 mL of Labeling Buffer (Component B).

Note 1: The unused Red Fluorescent Phalloidin Conjugate stock solution (Component A) should be aliquoted and stored at -20°C. Protect from light.

Note 2: Different cell types might be stained differently. The concentration of Red Fluorescent Phalloidin Conjugate working solution should be prepared accordingly.

B. Staining the Cells

Perform formaldehyde fixation. Incubate the cells with 3.0 –
 4.0 % formaldehyde in PBS at room temperature for 10 – 30 minutes.

Note: Avoid any methanol containing fixatives since methanol can disrupt actin during the fixation process. The preferred fixative is methanol-free formaldehyde

- 2. Rinse the fixed cells 2 3 times in PBS.
- 3. Optional: Add 0.1% Triton X-100 in PBS into fixed cells (from Step B.2) for 3 to 5 minutes to increase permeability. Rinse the cells 2–3 times in PBS.
- Add 100 μL/well (96-well plate) of 1x Red Fluorescent Phalloidin Conjugate working solution (from Step A) into the fixed cells (from Step B.2 or B.3), and stain the cells at room temperature for 15 – 60 minutes
- 5. Rinse cells gently with PBS 2 3 times to remove excess dye before plate sealing and imaging by using Texas Red channel.

ANTIBODY AND F-ACTIN COMBINATION STAINING

Note: This protocol is for one 96 - well plate.

A. Preparation of 1X Red Fluorescent Phalloidin Conjugate Solution

Add 10 µL Red Fluorescent Phalloidin Conjugate (Component A) to 10 mL of Labeling Buffer (Component B).

Note 1: The unused Red Fluorescent Phalloidin Conjugate solution (Component A) should be aliquoted and stored protected from light at -20 °C.

Note 2: Different cell types might be stained differently. The concentration of Red Fluorescent Phalloidin Conjugate working solution should be prepared accordingly.

B. Staining the Cells

Perform formaldehyde fixation. Incubate the cells with 3.0 –
 4.0 % formaldehyde in PBS at room temperature for 10–30 minutes.

Note: Avoid any methanol containing fixatives since methanol can disrupt actin during the fixation process. The preferred fixative is methanol-free formaldehyde

Rinse the fixed cells 2–3 times in PBS.

- 3. Permeabilize fixed cells with 0.2% Triton X-100 in PBS; incubate for 10 30 minutes to increase permeability. Rinse the cells once with PBS.
- 4. Add 100 μ I/well 5% Goat serum and leave for 30 minutes. Aspirate serum but do not rinse.
- 5. Add 100µl of primary antibody solution with 0.5% BSA in PBS and incubate for 1 hour. Make sure the entire are is covered.
- Rinse cells 3 times in PBS, each time for 5 minutes. Aspirate PBS.
- 7. Add Goat serum, incubate for 1 2 minutes and aspirate.
- 8. Add 100 μ L/well of secondary antibody solution diluted in 0.5% BSA in PBS buffer and 100 μ L/well 1X Red Fluorescent Phalloidin Conjugate (from Step A) working solution and stain the cells at room temperature for 30 60 minutes. Keep in the dark.
- If desired, add nuclear dye at relevant dilution in 0.5% BSA in PBS and incubate 10 – 30 minutes. Wash cells 2 – 3 times for 5 minutes in PBS.

Note: do not use PI as nuclear dye with this product as it emits in a very similar range and the signal might be masked.

- We recommend using DAPI as it emits on the blue or DRAQ5 (ab108410) as it emits on the far red.
- 10. Seal plate and examine samples in fluorescence microscope using red (Texas Red) channel (Ex/Em = 594/610 nm) for Factin staining. Use the appropriate channels to detect your secondary antibody and nuclear staining.

6. Data Analysis

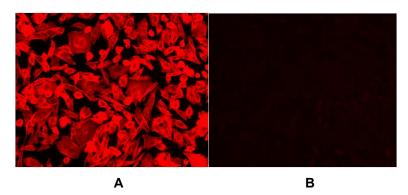


Figure 1. Image of CPA cells fixed with formaldehyde and stained with ab112126 in a black 96-well plate.

A: Cells were labeled with 1X Red Fluorescent Phalloidin Conjugate for 30 minutes only.

B: Cells were pre-treated with phalloidin for 10 minutes, then stained with 1X Red Fluorescent Phalloidin Conjugate for 30 minutes

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select "contact us" on www.abcam.com for the phone number for your region).



UK, EU and ROW

Email: technical@abcam.com | Tel: +44-

(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com | Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel: 01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 400 921 0189 / +86 21 2070 0500

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp

Copyright © 2013 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.