

Rhodamine Phalloidin Reagent ab235138

23 References

製品の概要

製品名	Rhodamine Phalloidin Reagent
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based (qualitative)
製品の概要	<p>Rhodamine Phalloidin Reagent ab235138 is one of a series of phalloidin conjugates that are used to stain actin filaments in fluorescence microscopy. The tetramethyl rhodamine dye can be easily detected with a fluorescent microscope at Ex/Em = 546/575 nm.</p> <p>Phalloidin conjugates are convenient probes for labeling, identifying and quantitating animal or plant actin filaments in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments.</p> <p>Review other popular phalloidin dye conjugates, including Phalloidin-iFluor 488, Phalloidin-iFluor 647, Phalloidin-iFluor 594, and Phalloidin-iFluor 555, search the website to see all phalloidin conjugates, or read the phalloidin staining protocol.</p>
特記事項	<p>Staining fixed cell or tissue samples with phalloidin conjugates is very simple; it requires a single 20-90 min incubation with the phalloidin, followed by 3 short wash steps. Phalloidin staining can be combined with antibody-based staining by adding the phalloidin conjugate during either the primary or secondary antibody incubation step.</p> <p>When used in unfixed samples, phalloidin binding leads to a decrease in the disassociation rate of actin subunits from the ends of actin filaments, essentially stabilizing actin filaments through the prevention of filament depolymerisation.</p>
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	300 tests
Rhodamine Phalloidin Conjugate	1 x 300 tests

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