

Nucleolar/Nuclear Staining Kit - Green/ Red Fluorescence - Cytopainter ab139474

[1 References](#) [画像数 1](#)

製品の概要

製品名	Nucleolar/Nuclear Staining Kit - Green/ Red Fluorescence - Cytopainter
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based
全工程の試験時間	0h 30m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Nucleolar/Nuclear Staining Kit - Green/ Red Fluorescence Cytopainter (ab139474) contains two proprietary dyes suitable for simultaneous live-cell staining of nucleoli and nuclei. The dyes allow examination of nucleolar dynamic changes in intracellular distribution, trafficking and localization arising from biological processes such as the cell cycle and ribosome biogenesis. The kit is compatible with most fluorescence detection systems, including conventional and confocal fluorescence microscopes, as well as High Content Screening (HCS) platforms.</p> <p>This kit is specifically designed for visualizing nucleoli and nuclei in living cells. A control nucleolus perturbation agent, actinomycin D, is provided for monitoring changes in nucleolar dynamics. Potential applications for this kit include monitoring impaired ribosome biogenesis, inhibition transcription, cell cycle dynamics and cellular stress, as well as the distribution, trafficking and dynamics of nucleolar proteins, the distribution of viral proteins, and potentially as an aid in identifying cancer cells.</p> <p>Review other dyes and kits for nucleolar staining, or the live cell staining fluorescent dyes guide</p>
試験プラットフォーム	Fluorescence microscope

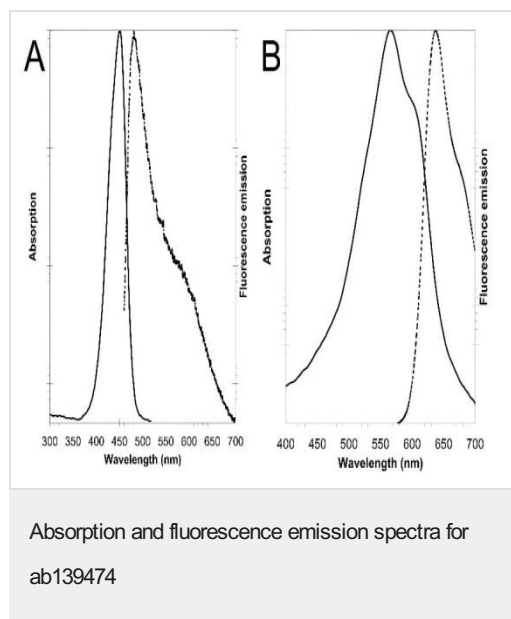
製品の特性

保存方法 Store at -80°C. Please refer to protocols.

内容	500 tests
10X Assay Buffer	1 x 15ml

内容	500 tests
Actinomycin D Control	1 x 125µg
Green Detection Reagent	1 x 50µl
Red Detection Reagent	1 x 50µl

画像



Absorption and fluorescence emission spectra for ab139474 Green (A) and Red (B) dyes. All spectra were determined in 1X Assay Buffer.

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