

# Nucleolar Staining Kit - Green Fluorescence - Cytopainter ab139475

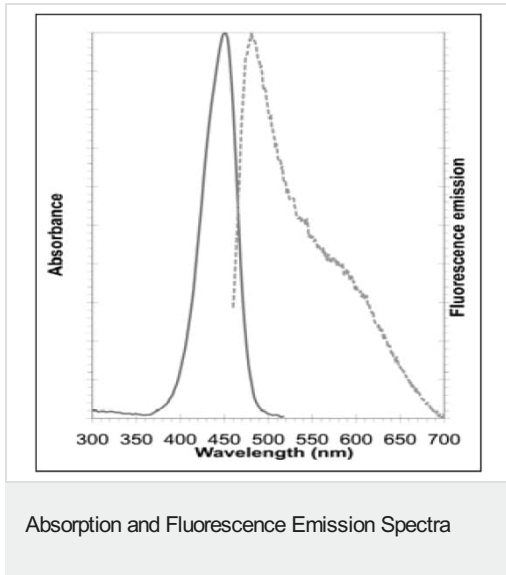
[1 Abreviews](#) [2 References](#) [1 图像](#)

### 概述

产品名称	Nucleolar Staining试剂盒- Green Fluorescence - Cytopainter
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based
检测时间	0h 30m
种属反应性	<b>与反应:</b> Mammals, Other species
产品概述	<p>Nucleolar Staining Kit - Green Fluorescence   Cytopainter (ab139475) contains a proprietary dye suitable for live-cell staining of nucleoli. The dye allows 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>ab139475 is specifically designed for visualizing nucleoli in living cells. The dye in the kit is resistant to photobleaching, facilitating its use in imaging applications. This nucleolar dye may be used in conjunction with common live-cell nuclear counterstains such as Hoechst 33342, Draq5™ or Nuclear Red dye.</p> <p>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 of 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 <a href="#">nucleolar staining</a>, or the <a href="#">live cell staining fluorescent dyes guide</a></p>
平台	Fluorescence microscope
性能	
存放说明	Please refer to protocols.

组件	500 tests
10X Assay Buffer	1 x 15ml
Actinomycin D Control	1 x 125µg
Green Detection Reagent	1 x 50µl

## 图片



Absorption and fluorescence emission spectra for Green Detection Reagent dye. Spectrum was 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.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