

# ER Staining Kit - Red Fluorescence - Cytopainter ab139482

★★★★☆ [4 Abreviews](#) [10 References](#) [画像数 2](#)

### 製品の概要

製品名	ER Staining Kit - Red Fluorescence - Cytopainter
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based
種交差性	<b>交差種:</b> Mammals, Other species
製品の概要	<p>ER Staining Kit - Red Fluorescence   Cytopainter (ab139482) contains an endoplasmic reticulum-selective red dye suitable for live cell, or detergent-permeabilized aldehyde-fixed cell staining. Micromolar concentrations of the red dye are sufficient for staining mammalian cells, as validated with human cervical carcinoma cell line HeLa, human T-lymphocyte cell line Jurkat and human bone osteosarcoma epithelial cell line, U2OS. (Ex/Em = 580/677 nm).</p> <p>One important application of the ER staining kit is in fluorescence co-localization imaging with green fluorescent protein (GFP)-tagged proteins, a powerful approach for determining the targeting of molecules to intracellular compartments and for screening of their associations and interactions.</p> <p>The ER staining kit is specifically designed for use with GFP-expressing cell lines, as well as cells expressing blue, cyan or yellow fluorescent proteins (BFPs, CFPs, YFPs). Additionally, the kit is suitable for use with live or post-fixed cells in conjunction with probes, such as labeled antibodies, or other fluorescent conjugates displaying similar spectral properties as fluorescein, or coumarin.</p> <p>A nuclear counterstain is also included in the kit.</p> <p>Review other dyes and kits for <a href="#">ER staining</a>, or the <a href="#">live cell staining fluorescent dyes guide</a></p>
特記事項	<p>This kit can be readily used in combination with other common UV and visible light excitable organic fluorescent dyes and various fluorescent proteins in multi-color imaging and detection applications. The dye emits in the Texas Red region of the visible light spectrum, and is resistant to photo-bleaching, concentration quenching and photoconversion.</p> <p>Previously called CytoPainter ER Staining Kit -Red Fluorescence.</p>
試験プラットフォーム	Fluorescence microscope

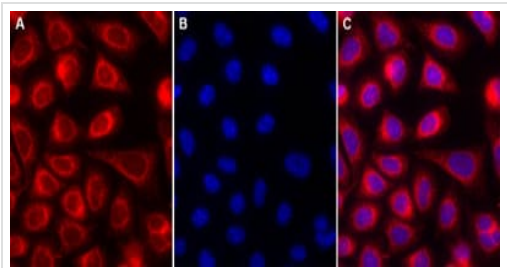
### 製品の特性

保存方法

Please refer to protocols.

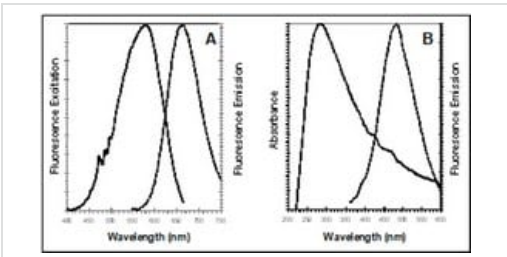
内容	500 tests
10X Assay Buffer	1 x 15ml
Hoechst 33342 Nuclear Stain	1 x 50µl
Red Detection Reagent	1 x 50µl

画像



Fluorescent staining using CytoPainter ER Staining kit

Live HeLa cells stained with Red Detection Reagent (A), Hoechst dye (B) and resulting composite image (C).



Fluorescence excitation and emission spectra for CytoPainter ER Staining kit

Fluorescence excitation and emission spectra for Red dye (panel A) and absorbance and fluorescent emission spectra for Hoechst 33342 dye (panel B). 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