

# Fixable Cell Viability Assay Kit (Fluorometric - Orange Ex 405 nm) - Cytopainter ab176740

画像数 2

## 製品概要

製品名	Fixable Cell Viability Assay Kit (Fluorometric - Orange Ex 405 nm) - Cytopainter
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 (quantitative)
種交差性	交差種: Mammals, Other species
製品の概要	Abcam's Fixable Cell Viability Assay Kit (Fluorometric - Orange Ex 405 nm)   Cytopainter (ab176740) is used to evaluate the viability of mammalian cells by flow cytometry.

This particular kit is designed to uniformly label fixed mammalian cells in blue fluorescence for flow cytometry applications with violet laser excitation. The kit uses a proprietary blue fluorescent dye that is more fluorescent upon bonding to cellular components. The kit provides all the essential components with an optimized cell-labeling protocol. It is an excellent tool for preserving of fluorescent images of particular cells, and can also be used for fluorescence flow cytometry applications.

The dye is designed to label cells at Ex/Em = 398/550 nm excited with a violet laser at 405 nm.

## 特記事項

### Related assays

Review the [cell health assay guide](#) to learn about kits to perform a [cell viability assay](#), [cytotoxicity assay](#) and [cell proliferation assay](#).

## 試験プラットフォーム

Flow cytometer

## 製品の特性

## 保存方法

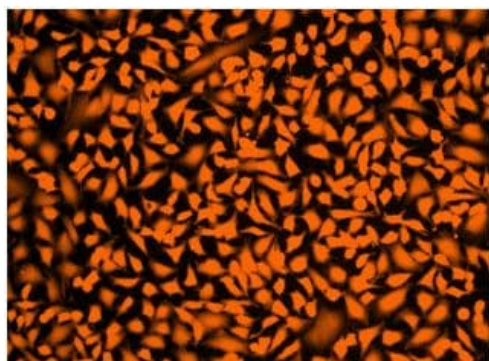
Store at -20°C. Please refer to protocols.

内容	200 tests
DMSO	1 x 200µl
Tracking dye Violet 550	1 vial

## 関連性

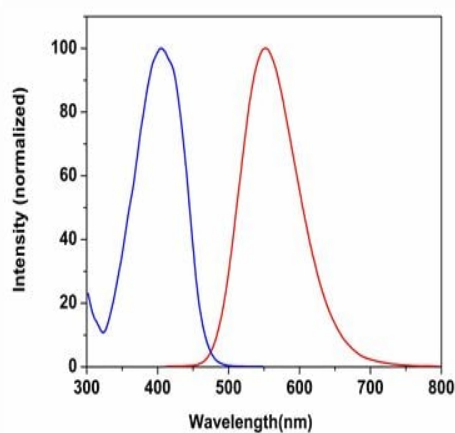
Cell viability is a determination of living or dead cells, based on a total cell population. Cell viability assess healthy cells in a sample, with no distinction between dividing or quiescent cells. An increase in cell viability indicates cell growth, while a decrease in viability can generally be interpreted as the result of either toxic effects of compounds/agents or suboptimal culture conditions.

## 画像



HeLa cells stained with CytoPainter Fixable Cell Viability Assay (Fluorometric-Orange Ex 405 nm) (ab176740)

HeLa cells fixed with formaldehyde and stained with Abcam's CytoPainter Fixable Cell Viability Assay Kit (Fluorometric - Orange Ex 405 nm) (ab176740) in a Costar black wall/clear bottom 96-well plate.



Excitation and Emission spectra for CytoPainter Fixable Cell Viability Assay Kit (Fluorometric - Orange Ex 405 nm) (ab176740)

Spectrum for ab176740.

**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