

ab176743

**Fixable Cell Viability Assay
Kit (Orange Fluorescence,
for 488 nm or 561 nm**

Instructions for Use

For evaluation of the viability of mammalian cells by fluorescent labeling and flow cytometry.

This product is for research use only and is not intended for diagnostic use.

Version: 3 Last Updated: 13 February 2026

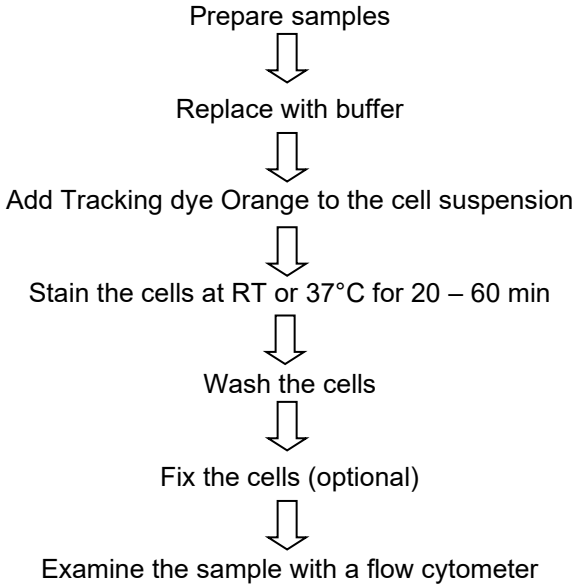
Table of Contents

Table of Contents	2
1. Introduction	3
2. Protocol Summary	4
3. Materials Supplied	4
4. Storage and Stability	5
5. Materials Required, Not Supplied	5
6. Assay Protocol	6
7. Data Analysis	8

1. Introduction

Abcam's Fixable Cell Viability Assay Kit (Orange Fluorescence, for 488 nm or 561 nm excitation) is used to evaluate the viability of mammalian cells by flow cytometry. The fluorescent dye provided in the kit is retained in cells by reacting with cellular components. For viable cells, only the cell-surface amines are available to react with the dye while for the necrotic cells or the other cells with compromised membranes, the reactive dye reacts with cell surface amines and intracellular amines, resulting in more intense fluorescent staining. The difference in fluorescence intensity between the live and dead cell populations is ~100-500 fold and can be completely preserved after fixation. The approximate fluorescence excitation is 547 nm and emission maximum is 573 nm. The Excitation source is 488 nm or 561 nm.

2. Protocol Summary



3. Materials Supplied

Item	200 tests
Tracking dye Orange	1 vial
DMSO	1 vial (200 µL)

4. Storage and Stability

Upon receipt, store kit at -20°C. Avoid exposure to light and moisture (store desiccated).

5. Materials Required, Not Supplied

- Sodium azide-free and serum/protein free buffer such as HHBS Buffer (1X Hanks and 20 mM HEPES buffer)
- CO₂ incubator
- Pipettes and pipette tips
- FACS tubes

6. Assay Protocol

1. Reagent Preparation:

a) 500X DMSO stock solution

Add 200 μ L DMSO to the vial of Tracking dye Orange.

NOTE: The unused stock solution should be aliquoted and stored at -20°C. Avoid repeated freeze/thaw cycles.

2. Sample Analysis:

a) Prepare cells for flow cytometry using 1X Hanks and 20 mM HEPES buffer (HHBS) or a sodium azide-free and serum/protein free buffer of your choice.

b) Wash cells once with HHBS or sodium azide-free and serum/protein free buffer of your choice.

c) Resuspend cells at 5-10 x 10⁶/mL in HHBS or sodium azide-free and serum/protein free buffer of your choice.

d) Add 1 μ L of Tracking dye Orange stock solution (see Reagent Preparation) to 0.5 mL of cells /assay and mix it well.

e) Incubate for 20-60 min at room temperature or 37°C, 5% CO₂ incubator, protected from light.

NOTE: The optimal stain concentrations and incubation time should be experimentally determined for different cell lines.

f) Wash cells twice and resuspend cells in with HHBS or buffer of your choice.

- g)** Fix cells as desired (optional).
- h)** Analyze cells with a flow cytometer.

7. Data Analysis

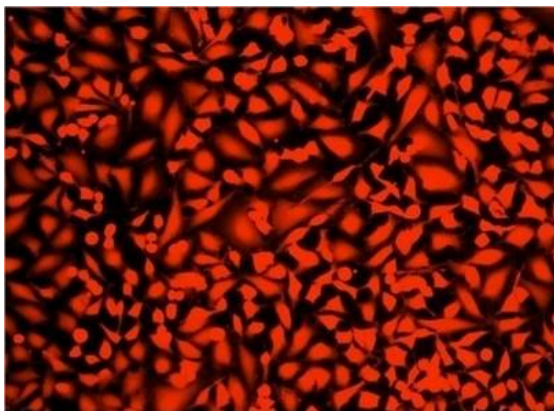


Figure 1. Fluorescent imaging of HeLa cells fixed with formaldehyde and labeled with ab176743 in a black wall/ clear bottom 96 well plate

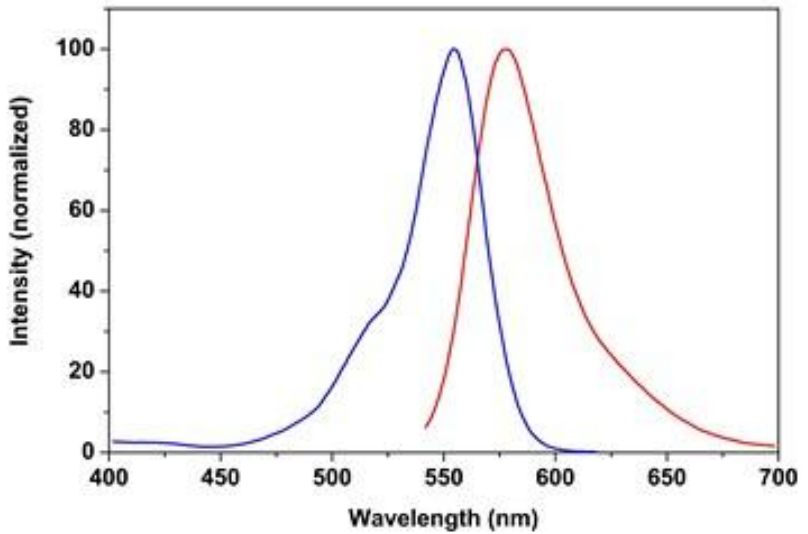


Figure 2. Excitation and Emission Spectra of Fixable Cell Viability Assay Kit (Orange Fluorescence, for 488 nm or 561 nm excitation) (ab176743).

Technical Support

Copyright © 2026 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)