

ab287858 – Live/Dead Cell Viability Assay Kit

For screening/studying/characterization of stimulators/inhibitors that affect cell viability.

For research use only - not intended for diagnostic use.

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab287858>

Storage and Stability

On receipt entire assay kit should be stored at -20°C, protected from light. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Materials Supplied

Item	Quantity	Storage Condition
Assay Buffer 27	100 mL	-20°C
Cell Dye II	1 vial	-20°C
Dead Cell Staining Dye	50 µL	-20°C

PLEASE NOTE: Assay Buffer 27 was previously labelled as Assay Buffer XXVII and Assay Buffer. The composition has not changed.

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- 6-well, 12-well, or 24-well plate
- 37°C Incubator with 5% CO₂
- Light and fluorescence microscope with Ex/Em = 485-495/530-635 nm
- FACS with Red and Green Channel detector

Reagent Preparation

- Read the entire protocol before performing the assay.
- Briefly spin the small vials prior to use.
- Open all reagents under sterile conditions (e.g. cell culture hood).

Assay Buffer 27: Store at -20°C. Warm to 37°C before use.

Cell Dye II: Reconstitute in 100 µL DMSO. Light sensitive, do not expose to intense light. Store at -20°C.

Dead Cell Staining Dye: Light sensitive, do not expose to intense light. Store at -20°C.

Assay Protocol

Δ Note: This protocol is for a 24-well plate. Adjust the volume according to the plate size.

Cell Culture and Staining:

1. Grow cells in 37°C incubator containing 5% CO₂ in desired media. Treat cells with compounds of interest, if desired. As a control, we recommend treating cells with vehicle alone.

Δ Notes:

- a) Adherent cells can be grown on cover slip for microscopy application to obtain better image resolution.

b) We recommend using suspension cells for flow cytometry application.

2. Mix 2 µL of Cell Dye II and 1 µL of Dead Cell Staining Dye in 1 mL of Assay Buffer 27. Prepare enough Staining Solution for your assay (0.5 mL per well in 24 well dish). Scale up accordingly for larger numbers of assays.
3. For suspension cells, collect ~1 x 10⁶ cells by centrifugation at 500 X g for 5 min. Resuspend in 0.5 mL Staining Solution. For adherent cells, remove the media carefully and add 0.5 mL Staining Solution to each well. Incubate for 15 min. at 37°C.

Detection:

Microscopy:

1. Place the cell suspension on a glass slide. Cover the cells with a glass coverslip.
2. For analyzing adherent cells, cell culture plates can be used directly. If cells are grown on a coverslip, invert coverslip on a glass slide and visualize cells.
3. Observe cells immediately under a light and fluorescence microscope (detects green and red wavelength [Ex/Em = 485-495/530-635 nm]). Cell Dye II stains healthy cells green. Dead Cell Staining Dye stains dead cell red.
4. Acquire several images per well for analysis.

Flow Cytometry:

1. Wash cells once with PBS.
2. Resuspend cell pellet in Assay Buffer 27 (~10⁶ cells/mL).
3. Analyze immediately using flow cytometry. Cell Dye II is measured in the FL1 channel and Dead Cell Staining Dye is measured in the FL3 channel.
4. To ensure that only proper target cells are gated, use a side scatter versus FL-1 plot.

Δ Notes:

- a) We recommend staining cells with Cell Dye II alone and Dead Cell Staining Dye alone to choose the proper instrument gating set up.
- b) We recommend keeping unstained control cells (i.e. without any Dye staining) suspended in Assay Buffer 27 for both treated and untreated samples to set up the flow cytometer instrument.

Technical Support

Copyright © 2025 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:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)