



ELISA PRODUCT INFORMATION & MANUAL

Live-Dead Cell Staining Kit (Fluorometric) (Fluorometric) *NBP2-54864*

Enzyme-linked Immunosorbent Assay for quantitative
detection. For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

Table of Contents

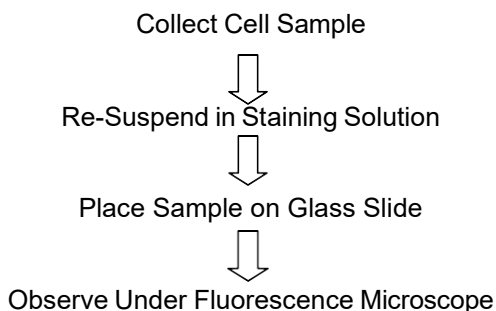
1.	Overview	3
2.	Protocol Summary	3
3.	Components and Storage	4
4.	Assay Protocol	5

1. Overview

Distinguishing between live and dead cells is very important for investigation of growth control and cell death. Novus Biologicals's Live-Dead Cell Staining Kit (Fluorometric) provides the ready-to-use reagents for convenient discrimination between live and dead cells.

The kit utilizes a cell-permeable green fluorescent dye (Ex/Em = 488/518 nm), to stain live cells. Dead cells can be easily stained by propidium iodide (PI), a cell non-permeable red fluorescent dye (Ex/Em = 488/615). Stained live and dead cells can be visualized by fluorescence microscopy using a band-pass filter (detects FITC and rhodamine). The kit provides sufficient reagents for 100 stainings using 24-well plate.

2. Protocol Summary



3. Components and Storage

A. Kit Components

Item	Quantity
Solution A (1 mM Live-Dye)	50 μ l
Solution B (2.5 mg/ml PI)	50 μ l
Staining Buffer	50 mL

* Store kit at -20°C. Protect from light. Store Staining Buffer at +4°C after opening. All reagents are stable for 1 year under proper storage conditions. Please read the entire protocol before performing the assay.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Fluorescence microscope
- Glass slides and coverslips
- Orbital shaker

4. Assay Protocol

1. Prepare enough Staining Solution for your assay (0.5 ml per well in 24 well dish):
Mix 1 μ l of Solution A and 1 μ l of Solution B in 1 ml of Staining Buffer. Scale up accordingly for larger numbers of assays.
2. Collect cells (1×10^6 cells) by centrifugation at $500 \times g$ for 5 min.
3. Re-suspend to 0.5 ml Staining Solution
4. Incubate for 15 min at 37°C .
5. Place the cell suspension on a glass slide. Cover the cells with a glass coverslip.

Note:

For analyzing **adherent cells**, grow cells directly on a coverslip. Following incubation with the Staining Solution, invert coverslip on a glass slide and visualize cells.

6. Observe cells immediately under a fluorescence microscope using a band-pass filter (detects fluorescein and rhodamine).

Healthy cells stain only the cell-permeable Live-Dye, fluorescing green. Dead cells can stain both the cell-permeable Live-Dye and the cell non-permeable PI (red), the overlay of green and red appears to be yellow-red.

Notes:

- a) As the optimal staining conditions may vary among different cell types, we recommend that a suitable concentration of Solution A and B be determined individually.
- b) Please note that PI is suspected to be highly carcinogenic, so careful handling of the reagent is required.