

# Live-Dead Cell Staining Kit

(Catalog #K501-100; 100 stainings; Store kit at -20°C)

## I. Introduction:

Distinguishing between live and dead cells is very important for investigation of growth control and cell death. The *Live-Dead Cell Staining Kit* provides the ready-to-use reagents for convenient discrimination between live and dead cells. The kit utilizes Live-Dye™, 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.

## II. Kit Contents:

Component	K501-100	Part Number
Solution A (1 mM Live-Dye)	50 µl	K501-100-1
Solution B (1 mg/ml PI)	50 µl	K501-100-2
Staining Buffer	50 ml	K501-100-3

## III. Cell Staining Protocol:

1. Prepare enough Staining Solution for your assay (0.5 ml per well in 24 well dish): mix 1 µl of Solution A and 1 µl of Solution B in 1 ml of Staining Buffer. Scale up accordingly for larger numbers of assays.
2. Collect cells (1 x 10<sup>6</sup> cells) by centrifugation at 500 X g for 5 min.
3. Resuspend 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. 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.

## IV. Caution:

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.

Please note that PI is suspected to be highly carcinogenic, so careful handling of the reagent is required.

## V. Storage and Stability:

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.

**FOR RESEARCH USE ONLY! Not to be used on humans.**

## RELATED PRODUCTS:

### Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents
- Apoptosis Inducers and Set
- Apoptosis siRNA Vectors

### Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Rapid Separation Kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

### Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- High Throughput Apoptosis/Cell Viability Assay Kits
- LDH-Cytotoxicity Assay Kit
- Bioluminescence Cytotoxicity Assay Kit
- Live/Dead Cell Staining Kit

### Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

### Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits
- Beta-Secretase Activity Assay Kit

### Adipocyte & Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP Activity Assay & Drug Discovery Kits
- PLTP Activity Assay & Drug Discovery Kits
- Total Cholesterol Quantification Kit

### Molecular Biology & Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Min DNA Ligation Kit
- 20 Min Gel Staining/Destaining Kit
- β-Galactosidase Staining Kit & Luciferase Reporter Assay Kit

### Growth Factors and Cytokines

### Monoclonal and Polyclonal Antibodies