

Live and Dead Cell Assay ab115347

[47 References](#) [4 Images](#)

Overview

Product name	Live and Dead Cell Assay
Detection method	Fluorescent
Product overview	<p>Live Dead Assay Kit ab115347 differentially labels live and dead cells with fluorescent dyes with a one-step live dead assay protocol. It is used for the rapid quantitation of cell viability using flow cytometry or fluorescent microscopy.</p> <p>The Live Dead assay staining solution is a mixture of two fluorescent dyes that differentially label live and dead cells.</p> <p>The Live cell dye labels intact, viable cells green. It is membrane permeant and non-fluorescent until ubiquitous intracellular esterases remove ester groups and render the molecule fluorescent. The Excitation (max) and Emission(max) are 494nm and 515nm (similar to FITC).</p> <p>The Dead cell dye labels cells with compromised plasma membranes red. It is membrane-impermeant and binds to DNA with high affinity. Once bound to DNA, the fluorescence increases >30-fold. The Excitation (max) and Emission(max) are 528nm and 617nm.</p> <p>The Live Dead assay protocol uses a one-step staining procedure that is simple and fast. It can be used directly in cell culture media.</p>
Notes	<p>This assay is not suitable for use with fixed cells / cell fixation.</p> <p>The Live Dead assay staining solution provided is sufficient for ~1000 assays.</p> <p>Related assays</p> <p>Review the cell health assay guide to learn about kits to perform a cell viability assay, cytotoxicity assay and cell proliferation assay.</p>
Platform	Flow cytometer, Fluorescence microscope

Properties

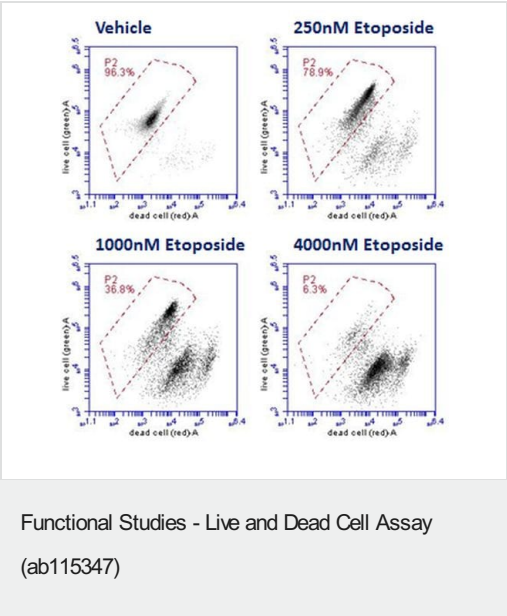
Storage instructions Store at -20°C. Please refer to protocols.

Components	1000 tests
1000X Live/Dead Cell stain in DMSO	1 x 0.1ml

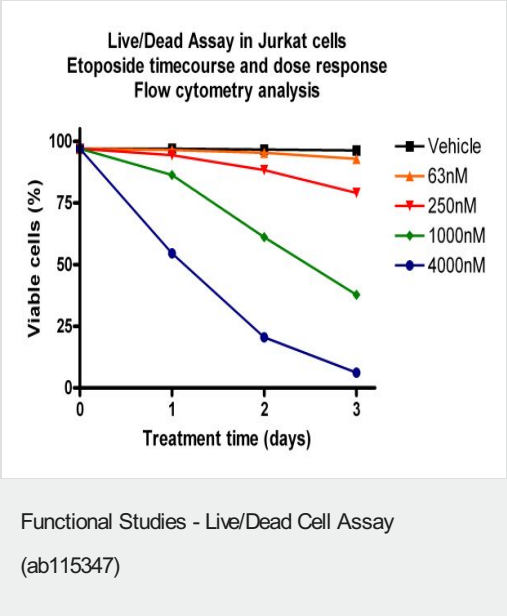
Relevance

Distinguishing between live and dead cells is very important for investigation of growth control and cell death.

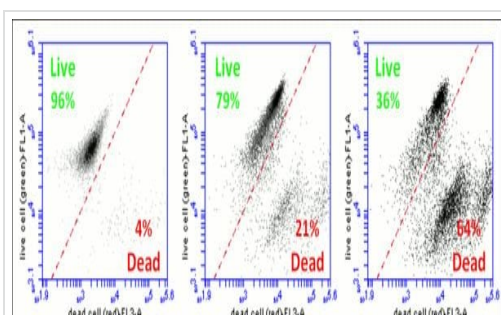
Images



Dot plots showing live/dead analysis of vehicle or drug treated Jurkat cells (day 3 of treatment). The indicated drug is used to induce cell death. Live cells are on the y-axis and dead cells are on the x-axis. The red polyongate identifies live cells and the number indicates the percent of live cells.



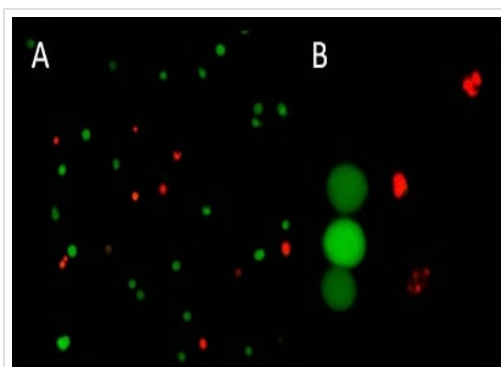
Quantification of % viable cells of Jurkat cells treated with a dose response of the inidicated drug (to induce cell death) and analyzed using the live/dead assay stain on days 1, 2 and 3 using flow cytometry.



Functional Studies - Live/Dead Cell Assay

(ab115347)

The sample dot plots demonstrate varying ratios of live and dead cells. More green = upper left = live cells; more red = lower right = dead cells.



Functional Studies - Live/Dead Cell Assay

(ab115347)

Jurkat cells stained with the live/dead assay kit. Jurkat cells treated with a drug to induce cell death were labeled with the live/dead assay stain. Live cells (with esterase activity) stain green and dead cells (compromised plasma membrane) stain red. (A) Field of cells following 10 minute staining in media of live/dead stain. (B) Magnified view showing that in live cells the whole cell is stained green whereas in dead red cells it is the fragmented nuclear DNA that is stained.

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors