

Apoptosis and Necrosis Quantitation Kit Plus

A convenient assay for quantifying apoptotic (green) and necrotic (red) cells within the same cell population by flow cytometry or fluorescence microscopy.

Biodum

Product Description

The Apoptosis and Necrosis Quantitation Kit Plus provides a convenient assay for quantifying apoptotic (green) and necrotic (red) cells within the same cell population by flow cytometry or fluorescence microscopy.

Apoptosis and necrosis are two major processes by which cells die. Apoptosis is an active, genetically regulated disassembly of the cell from within. During apoptosis, phosphatidylserine (PS) is translocated from the inner to the outer leaflet of the plasma membrane for phagocytic cell recognition. The human anticoagulant Annexin V is a 35 kDa, Ca²⁺-dependent phospholipid-binding protein with high affinity for PS. Annexin V labeled with CF®488A stains apoptotic cells green by binding to PS exposed on the cell surface. CF®488A is spectrally similar to fluorescein (FITC), with much brighter and more photostable fluorescence.

Necrosis normally results from severe cellular insult. Both internal organelle and plasma membrane integrity are lost, resulting in the release of cellular contents into the surrounding environment. Ethidium Homodimer III (EthD-III) is a highly positively charged nucleic acid probe, which is impermeant to live cells, but stains cells with compromised membrane integrity with red fluorescence. EthD-III is a superior alternative to Propidium Iodide (PI) or Ethidium Homodimer I because of its significantly higher affinity for DNA and higher fluorescence quantum yield. The Apoptosis & Necrosis Quantitation Kit Plus provides a convenient assay for quantifying apoptotic (green) and necrotic (red) cells within the same cell population by flow cytometry or fluorescence microscopy.

See our full selection of <u>Cell Viability & Apoptosis Assays</u> and <u>Viability & Apoptosis Assays for Flow Cytometry</u>.

CF dye technology is covered by US and international patents.

This datasheet was generated on November 20, 2025 at 09:35:47 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/apoptosis-and-necrosis-quantitation-kit-plus/

Product attributes

Call us: 800-304-5357

Apoptosis/viability marker	Phosphatidylserine/Annexin V, Dead cell stain, Apoptosis/necrosis assay
For live or fixed cells	For live/intact cells
Detection method/readout	Fluorescence microscopy, Flow cytometry
Assay type/options	Endpoint assay
Colors	Green/Red
Product origin	Annexin V (human); recombinant, produced in E. coli
Storage Conditions	Store at 2 to 8 °C, Do not freeze, Protect from light