

## Product datasheet

# Apoptosis/ Necrosis Assay Kit (blue, red, green) ab176750

★★★★★ [1 Abreviews](#) [13 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Apoptosis/ Necrosis Assay Kit (blue, red, green)
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based
<b>Assay time</b>	1h 00m
<b>Product overview</b>	Apoptosis/ Necrosis Assay Kit (blue, red, green) (ab176750) is designed to simultaneously monitor apoptotic, necrotic and healthy cells.

The PS sensor used in this kit has red fluorescence (Ex/Em = 630/660 nm) upon binding to membrane PS.

Necrosis has been characterized as passive, accidental cell death resulting from environmental perturbations with uncontrolled release of inflammatory cellular contents.

Loss of plasma membrane integrity, as demonstrated by the ability of the membrane-impermeable DNA Nuclear Green DCS1 dye (Ex/Em = 490/525 nm) to label the nucleus, represents a straightforward approach to demonstrate late stage apoptosis and necrosis.

This apoptosis / necrosis assay also provides a live cell cytoplasm labeling dye CytoCalcein Violet 450 (Ex/Em = 405/450 nm) to label live cell cytoplasm.

This kit is optimized to simultaneously detect cell apoptosis (Red), apoptosis (red and/or green) and healthy cells (blue) with a flow cytometer or fluorescence microscope.

### Notes

This product was previously called Apoptosis/ Necrosis Detection Kit.

### Other apoptosis assays

For more apoptosis assays, review the [apoptosis assay and apoptosis marker guide](#).

### Platform

Flow cytometer, Fluorescence microscope

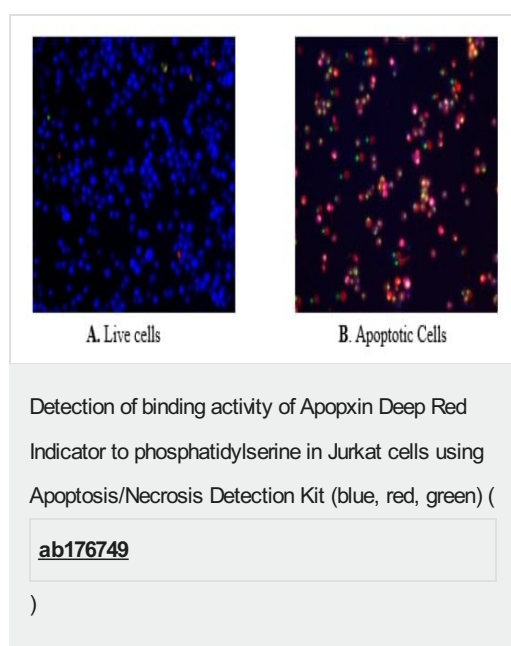
### Properties

## Storage instructions

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

Components	100 tests
Apopxin Deep Red Indicator	1 x 200µl
Assay Buffer	1 x 50ml
CytoCalcein Violet 450	1 vial
Nuclear Green DCS1 Dye	1 x 100µl

## Images



The fluorescence image shows cells that are live (blue, stained by CytoCalcein Violet 450), apoptotic (red, stained by Apopxin Deep Red Indicator), and necrotic (green, indicated by Nuclear Green DCS1 staining) in Jurkat cells induced by 1µM staurosporine for 3 hours. The fluorescence images of the cells were taken with a fluorescence microscope through the Violet, Cy5 and FITC channel respectively. Individual images taken from each channel from the same cell population were merged as shown above. A: Non-induced control cells; B: Triple staining of staurosporine-induced cells.

**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