



Caspase-3 Staining Kit (Red)

Catalog Number KA0744

96 assays

Version: 03

Intended for research use only

Introduction and Background

A. Overview

Activation of caspases plays a central role in apoptosis. The Red Active Caspase Staining Kit provides a convenient means for detecting activated caspases-3 in living cells. The assay utilizes a caspase-3 inhibitor DEVD-FMK conjugated to sulfo-rhodamine (Red-DEVD-FMK) as the fluorescent in situ marker. Red-DEVD-FMK is cell permeable, nontoxic, and irreversibly binds to activated caspases-3 in apoptotic cells. The red fluorescence label allows for direct detection of activated caspases-3 in apoptotic cells by fluorescence microscopy, flow cytometry, or fluorescence plate reader.

Material and Method

A. List of component

1. Red-DEVD-FMK 100 μ l.
2. Wash Buffer 2 x 100 ml.
3. Z-VAD-FMK 10 μ l.

B. Stability and storage

Store kit at -20°C.

C. Protocol

1. Staining:
 - a. Induce apoptosis in cells (1×10^6 /ml) by desired method. Concurrently incubate a control culture without induction. An additional control can be prepared by adding the caspase inhibitor Z-VAD-FMK at 1 μ l/ml to an induced culture to inhibit caspase activation.
 - b. Aliquot 300 μ l each of the induced and control cultures into eppendorf tubes.
 - c. Add 1 μ l of Red-DEVD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO₂.
 - d. Centrifuge cells at 3000 rpm for 5 minutes and remove supernatant.
 - e. Resuspend cells in 0.5 ml of Wash Buffer, and centrifuge again.
 - f. Repeat Step e.
 - g. Proceed to 2, 3, or 4 depending on methods of analysis.
2. Quantification by Flow Cytometry:

For flow cytometric analysis, resuspend cells in 300 μ l of Wash buffer. Put samples on ice. Analyzing samples by flow cytometry using the FL-2 channel.
3. Detection by Fluorescence Microscopy:

For fluorescence microscopic analysis, resuspend cells in 100 μ l Wash Buffer. Put one drop of the cell suspension onto a microslide and cover with a coverslip. Observe cells under a fluorescence microscope

using rhodamine filter. Caspase positive cells appear to have brighter green signals, whereas caspase negative control cells show much weaker signal.

4. Analysis by Fluorescence Plate Reader:

For analysis with fluorescence plate reader, resuspend cells in 100 μ l Wash Buffer and then transfer the cell suspension to each well of the black microtiter plate. Measure the fluorescence intensity at Ex/Em = 540/570 nm (Note: Ex/Em = 488/570 nm will also work, although it's not an optimal wavelength). For control, use wells containing unlabeled cells.