

Introduction and Background

A. Overview

Activation of caspases plays a central role in apoptosis. The Fluorescein Active Caspase-2 Staining Kit provides a convenient means for sensitive detection of activated caspase-2 in living cells. The assay utilizes the caspase-2 inhibitor, VDVAD-FMK, conjugated to FITC (FITC-VDVAD-FMK) as a marker. FITCVDVAD-FMK is cell permeable, nontoxic, and irreversibly binds to activated caspase-2 in apoptotic cells. The FITC label allows for direct detection of activated caspases in apoptotic cells by fluorescence microscopy, flow cytometry, or fluorescence plate reader.

B. Notice for Application of Kit

- ✓ This kit has been configured for research use only and is not for diagnostic and clinical use.

Material and Method

A. List of component

1. FITC-VDVAD-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 negative control can be prepared by adding the caspase inhibitor Z-VAD-FMK at 1 μ l/ml to an induced culture to inhibit caspase-2 activation.
 - b. Aliquot 300 μ l each of the induced and control cultures into eppendorf tubes.
 - c. Add 1 μ l of FITC-VDVAD-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. Keep samples on ice. Analyzing samples by flow cytometry using the FL-1 channel.

3. Detection by Fluorescence Microscopy:

For fluorescence microscopic analysis, resuspend cells in 100 μ l Wash Buffer. Transfer one drop of the cell suspension onto a microslide and cover with a coverslip. Observe cells under a fluorescence microscope using FITC 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 into each well in the black microtiter plate. Measure the fluorescence intensity at Ex. = 485 nm and Em. = 535 nm. For control, use wells containing unlabeled cells.