

Instruction Manual

Catalog Number	PK-CA577-K180-25	
Description	Activation of caspases plays a central role in apoptosis. The PromoKine Green Multi-Caspase Staining Kit provides a convenient means for sensitive detection of activated caspases in living cells. The assay utilizes the caspase family inhibitor, VAD-FMK, conjugated to FITC (FITC-VAD-FMK) as a marker. FITC-VAD-FMK is cell-permeable, nontoxic, and irreversibly binds to activated caspases 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.	
Quantity	25 assays	
Kit Components	Components	Quantity
	FITC-VAD-FMK	25 μ l
	Wash Buffer	50 ml
	Z-VAD-FMK	10 μ l
Applications / Assay Protocol	<p>A. Staining Procedure:</p> <ol style="list-style-type: none"> 1. 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 μl/ml to an induced culture to inhibit Caspase-3 activation. 2. Aliquot 300 μl each of the induced and control cultures into eppendorf tubes. 3. Add 1 μl of FITC-VAD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO₂. 4. Centrifuge cells at 3,000 rpm for 5 minutes and remove supernatant. 5. Resuspend cells in 0.5 ml of Wash Buffer, and centrifuge again. 6. Repeat Step 5. <p>Proceed to B, C, or D depending on methods of analysis.</p> <p>B. Quantification by Flow Cytometry: For flow cytometric analysis, resuspend cells in 300 μl of Wash buffer. Keep samples on ice. Analyze samples by flow cytometry using the FL-1 channel.</p> <p>C. Detection by Fluorescence Microscopy: For fluorescence microscopic analysis, resuspend cells in 100 μl Wash buffer. Put 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.</p> <p>D. Analysis by Fluorescence Plate Reader: For analysis with fluorescence plate reader, resuspend cells in 100 μl Wash Buffer and then transfer the cell suspension to 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.</p>	
Storage & Stability	Store kit at -20°C upon arrival. Store individual reagents as indicated on the respective labels.	

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.