

Instruction Manual

Catalog Number	PK-CA577-K199-100	
Description	Activation of caspases plays a central role in apoptosis. The PromoKine Red Caspase-9 Staining Kit provides a convenient means for detecting activated Caspase-9 in living cells. The assay utilizes a Caspase-9 inhibitor LEHD-FMK conjugated to sulfo-rhodamine (Red-LEHD-FMK) as the fluorescent marker. Red-LEHD-FMK is cell permeable, nontoxic, and irreversibly binds to activated Caspase-9 in apoptotic cells. The fluorescence label allows detection of activated Caspase-9 in apoptotic cells by fluorescence microscopy, flow cytometry, or using a fluorescence plate reader.	
Quantity	100 assays	
Kit Components	Components	Quantity
	Red-LEHD-FMK	100 µl
	Wash Buffer	2 x 100 ml
	Z-VAD-FMK	10 µl
Applications / Assay Protocol	<p>A. Staining Procedure</p> <ol style="list-style-type: none"> 1. Induce apoptosis in cells (1 x 10⁶/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-9 activation. 2. Aliquot 300 µl each of the induced and control cultures into eppendorf tubes. 3. Add 1 µl of Red-LEHD-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 <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. Analyzing samples by flow cytometry using the FL-2 channel.</p> <p>C. Detection by Fluorescence Microscopy: For fluorescence microscopic analysis, resuspend cells in 100 µl Wash Buffer. Transfer one drop of the cell suspension onto a microslide and cover with a coverslip. Observe cells under a fluorescence microscope using a rhodamine filter. Caspase-9 positive cells appear to have brighter green signals, whereas Caspase-9 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 into each well of a black microtiter plate. Measure the fluorescence intensity at Ex. = 540 nm and Em. = 570 nm (Note: Ex = 488 nm and Em. = 570 nm will also work, although it's not an optimal wavelength). 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.