

## Instruction Manual

<b>Catalog Number</b>	PK-CA577-K189-25									
<b>Description</b>	Activation of caspases plays a central role in apoptosis. The PromoKine Green Caspase-9 Staining Kit provides a convenient means for sensitive detection of activated Caspase-9 in living cells. The assay utilizes the Caspase-9 inhibitor LEHD-FMK conjugated to FITC (FITC-LEHD-FMK) as a fluorescent marker. FITC-LEHD-FMK is cell permeable, nontoxic, and irreversibly binds to activated Caspase-9 in apoptotic cells. The FITC label allows detection of activated caspases in apoptotic cells directly by fluorescence microscopy, flow cytometry, or using a fluorescence plate reader.									
<b>Quantity</b>	25 assays									
<b>Kit Components</b>	<table border="1"> <thead> <tr> <th>Components</th> <th>Quantity</th> </tr> </thead> <tbody> <tr> <td>FITC-LEHD-FMK</td> <td>25 µl</td> </tr> <tr> <td>Wash Buffer</td> <td>50 ml</td> </tr> <tr> <td>Z-VAD-FMK</td> <td>10 µl</td> </tr> </tbody> </table>	Components	Quantity	FITC-LEHD-FMK	25 µl	Wash Buffer	50 ml	Z-VAD-FMK	10 µl	
Components	Quantity									
FITC-LEHD-FMK	25 µl									
Wash Buffer	50 ml									
Z-VAD-FMK	10 µl									
<b>Applications / Assay Protocol</b>	<p><b>A. Staining Procedure</b></p> <ol style="list-style-type: none"> <li>1. Induce apoptosis in cells (<math>1 \times 10^6</math>/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.</li> <li>2. Aliquot 300 µl each of the induced and control cultures into eppendorf tubes.</li> <li>3. Add 1 µl of FITC-LEHD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO<sub>2</sub>.</li> <li>4. Centrifuge cells at 3,000 rpm for 5 minutes and remove supernatant.</li> <li>5. Resuspend cells in 0.5 ml of Wash Buffer, and centrifuge again.</li> <li>6. Repeat Step 5. Proceed to B, C, or D depending on methods of analysis.</li> </ol> <p><b>B. Quantification by Flow Cytometry:</b> For flow cytometric analysis, resuspend cells in 300 µl of Wash Buffer. Keep samples on ice. Analyzing samples by flow cytometry using the FL-1 channel.</p> <p><b>C. Detection by Fluorescence Microscopy:</b> 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 FITC filter. Caspase-9 positive cells appear to have brighter green signals, whereas Caspase-9 negative control cells show much weaker signal.</p> <p><b>D. Analysis by Fluorescence Plate Reader:</b> 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. = 485 nm and Em. = 535 nm. For control, use wells containing unlabeled cells.</p>									
<b>Storage &amp; Stability</b>	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.