



Caspase-8 Staining Kit (Red)

Catalog Number KA0760

100 assays

Version: 02

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Background	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Precaution of Use	4
Assay Protocol	5
Assay Procedure	5

Introduction

Background

Activation of caspases plays a central role in apoptosis. The Red Active Caspase-8 Staining Kit provides a convenient means for detecting activated caspase-8 in living cells. The assay utilizes a caspase-8 inhibitor IETD-FMK conjugated to sulfo-rhodamine (Red-IETD-FMK) as the fluorescent marker. Red-IETD-FMK is cell permeable, nontoxic, and irreversibly binds to activated caspase-8 in apoptotic cells. The fluorescence label allows detection of activated caspase-8 in apoptotic cells directly by fluorescence microscopy, flow cytometry, or fluorescence plate reader.

General Information

Materials Supplied

List of component

Component	Amount
Red-IETD-FMK	100 µl
Wash Buffer	2 x 100 ml
Z-VAD-FMK	10 µl

Storage Instruction

Store kit at -20 °C.

Precaution of Use

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

Assay Protocol

Assay Procedure

- Staining:
 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 μ l/ml to an induced culture to inhibit caspase activation.
 2. Aliquot 300 μ l each of the induced and control cultures into eppendorf tubes.
 3. Add 1 μ l of Red-IETD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO₂.
 4. Centrifuge cells at 3000 rpm for 5 minutes and remove supernatant.
 5. Resuspend cells in 0.5 ml of Wash Buffer, and centrifuge again.
 6. Repeat Step 5.
Proceed to 2, 3, or 4 depending on methods of analysis.
- 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.
- 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-8 positive cells appear to have brighter red signals, whereas caspase-8 negative control cells show much weaker signal.
- Analysis by Fluorescence Plate Reader:

For analysis with fluorescence plate reader, resuspend cells in 100 μ l Wash Buffer and 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.