



ELISA PRODUCT INFORMATION & MANUAL

Active Caspase Assay Kit (Fluorometric)

NBP2-54857

Enzyme-linked Immunosorbent Assay for quantitative
detection. For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

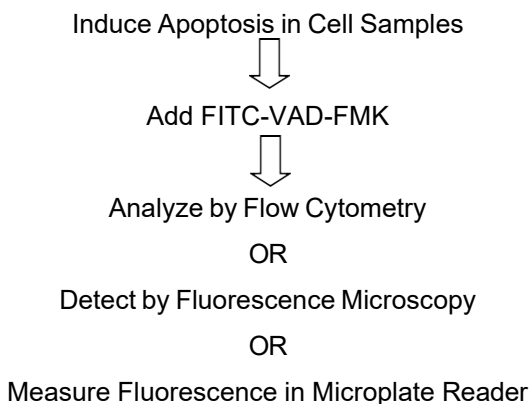
Table of Contents

1.	Overview	2
2.	Protocol Summary	2
3.	Components and Storage	3
4.	Assay Protocol	4
5.	Troubleshooting	6

1. Overview

Activation of Caspases plays a central role in apoptosis. Novus Biologicals's Active Caspase Assay Kit (Fluorometric) provides a convenient and sensitive means for detecting 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.

2. Protocol Summary



3. Components and Storage

A. Kit Components

Item	Quantity
FITC-VAD-FMK	100 μ L
Wash Buffer	2 x 100 mL
Z-VAD-FMK	10 μ L

* Store kit at -20°C.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Fluorescent microplate reader or microscope
- Flow cytometer
- Black microtiter plate
- Orbital shaker

4. Assay Protocol

1. Sample Preparation:

- a) Induce apoptosis in cells ($1 \times 10^6/\text{ml}$) 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 $1 \mu\text{l}/\text{ml}$ to an induced culture to inhibit Caspase activation.

Notes:

This product detects proteolytic activity. Do not use protease inhibitors in the sample preparation step as it might interfere with the assay.

- b) Aliquot $300 \mu\text{l}$ each of the induced and control cultures into eppendorf tubes.
- c) Add $1 \mu\text{l}$ of FITC-VAD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO_2 .
- d) Centrifuge cells at 3000 rpm for 5 minutes and remove supernatant.
- e) Re-suspend cells in 0.5 ml of Wash Buffer, and centrifuge again.
- f) Repeat Step e.

g) Proceed to Step 2, 3 or 4 depending on methods of analysis.

2. Quantification by Flow Cytometry:

For flow cytometric analysis, re-suspend cells in 300 μ l of Wash Buffer. Keep samples on ice. Analyze samples by flow cytometry using the FL-1 channel.

3. Detection by Fluorescence Microscopy:

For fluorescence microscopic analysis, re-suspend 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 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, re-suspend cells in 100 μ l Wash Buffer and then transfer the cell suspension to each well in the black microtiter plate.

Measure the fluorescence intensity at Excitation = 485 nm and Emission = 535 nm. For control, use wells containing unlabeled cells.

5. Troubleshooting

Problem	Reason	Solution
High Background	Cell density is higher than recommended	Refer to datasheet and use the suggested cell number
	Increased volumes of components added	Use calibrated pipettes accurately
	Incubation of cell samples for extended periods	Refer to datasheets and incubate for exact times
	Use of extremely confluent cells	Perform assay when cells are at 80-95% confluency
	Contaminated cells	Check for bacteria/ yeast/ mycoplasma contamination
Lower signal levels	Cells did not initiate apoptosis	Determine the time-point for initiation of apoptosis after induction (time-course experiment)
	Very few cells used for analysis	Refer to data sheet for appropriate cell number
	Incorrect setting of the equipment used to read samples	Refer to datasheet and use the recommended filter setting
	Use of expired kit or improperly stored reagents	Always check the expiry date and store the components appropriately

Erratic results	Uneven number of cells seeded in the wells	Seed only healthy cells (correct passage number)
	Adherent cells dislodged at the time of experiment	Perform experiment gently and in duplicates or triplicates for each treatment
	Incorrect incubation times or temperatures	Refer to datasheet & verify correct incubation times and temperatures
	Incorrect volumes used	Use calibrated pipettes and aliquot correctly

