

**ab65617**

# **Caspase-3 (active) Red Staining Kit**

## Instructions for Use

For the rapid, sensitive and accurate detection of activated Caspase-3 in living cells.

This product is for research use only and is not intended for diagnostic use.

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

# Table of Contents

---

1. Overview	2
2. Protocol Summary	2
3. Components and Storage	3
4. Assay Protocol	4
5. Factors to consider for caspase activity assays	6
6. Troubleshooting	8

## 1. Overview

---

Activation of caspases plays a central role in apoptosis. Abcam's Caspase-3 (active) Red Staining Kit provides a convenient means for detecting activated Caspase-3 in living cells. The assay utilizes the Caspase-3 inhibitor DEVD-FMK conjugated to sulfo-rhodamine (Red-DEVD-FMK) as the fluorescent *in situ* marker. Red-DEVD-FMK is cell permeable, nontoxic, and irreversibly binds to activated Caspase-3 in apoptotic cells.

## 2. Protocol Summary

---

Sample Preparation



Analyze by Flow Cytometry

OR

Detect by Fluorescence Microscopy

OR

Measure Fluorescence in Microplate Reader



### 3. Components and Storage

---

#### A. Kit Components

Item	Quantity
Red-DEVD-FMK	100 $\mu$ L
Wash Buffer IV/Wash Buffer	2 x 100 mL
Z-VAD-FMK	10 $\mu$ L

\* Store kit at  $-20^{\circ}\text{C}$ .

#### 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. Staining Procedure

- a) Induce apoptosis in cells ( $1 \times 10^6$  cells/ml) by the desired method. Concurrently incubate a control culture *without* induction.

An additional **control** can be prepared by adding the caspase inhibitor Z-VAD-FMK at 1  $\mu$ l/ml to an induced culture to inhibit caspase activation.

#### Note:

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  $\mu$ l each of the induced and control cultures into eppendorf tubes.
- c) Add 1  $\mu$ l of Red-DEVD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO<sub>2</sub>.
- d) Centrifuge cells at 3000 rpm for 5 minutes and remove supernatant.
- e) Re-suspend cells in 0.5 ml of Wash Buffer IV/Wash Buffer, and centrifuge again.

- f) Repeat step e.
- g) Proceed to Step 2, 3 or 4 depending on methods of analysis of the un-induced control.

### **2. Quantification by Flow Cytometry:**

For flow cytometric analysis, re-suspend cells in 300 µl of Wash Buffer IV/Wash buffer. Put samples on ice. Analyze samples by flow cytometry using the FL-2 channel.

### **3. Detection by Fluorescence Microscopy:**

For fluorescence microscopic analysis, re-suspend cells in 100 µl Wash Buffer IV/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 positive cells appear to have brighter red 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 µl Wash Buffer IV/Wash Buffer and then transfer the cell suspension to each well of the black microtiter plate. Measure the fluorescence intensity at Ex/Em = 540/570 nm.

For control, use wells containing unlabeled cells.

### **Note:**

Ex/Em = 488/570 nm will also work, although it's not an optimal wavelength.

## **5. Factors to consider for caspase activity assays**

---

Three major factors need to be taken into account when using caspase activity assays:

1. The substrate in a particular assay is not necessarily specific to a particular caspase.

Cleavage specificities overlap so reliance on a single substrate/assay is not recommended. Other assays, such as Western blot or use of fluorescent substrates e.g. FRET assays should be used in combination with caspase activity assays.

2. The expression and abundance of each caspase in a particular cell type and cell line will vary.
3. As the activation and cleavage of caspases in the cascade will change over time, you should consider when particular caspase will be at its peak concentration e.g. after 3 hours, after 20 hours etc.

The table below shows the known cross-reactivities with other caspases.

Classification of caspases based on synthetic substrate preference, does not reflect the real caspase substrate preference in vivo and may provide inaccurate information for discriminating amongst



caspase activities. Thus, caution is advised in applying the intrinsic tetrapeptide preferences to predict the targets of individual caspases.

### Apoptotic Executer Caspases

Caspase	Cleavage motif	Inhibitor motif	Cross-reactivity with other caspase:											
			1	2	3	4	5	6	7	8	9	10		
<b>Caspase -3</b>	DEVD	DEVD, LEHD*, IETD, LETD		Y						Y				
<b>Caspase 6</b>	VEID	DEVD, LEHD*, IETD, LETD			Y									
<b>Caspase 7</b>	DEVD	DEVD, LEHD*, IETD, LETD	Y		Y									

*\* inhibits at high concentration*

## 6. Troubleshooting

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
<b>High Background</b>	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
<b>Lower signal levels</b>	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

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
<b>Erratic results</b>	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



## **Technical Support**

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)