

ab65619

Caspase 9 (active) Red Staining Kit

Instructions for Use

For the rapid, sensitive and accurate detection of activated Caspase 9 in living cells

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

Table of Contents

| | |
|--|---|
| 1. Overview | 1 |
| 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 9 (active) Red Staining Kit provides a convenient means for detecting activated caspase 9 in living cells.

The assay utilizes the caspase 9 inhibitor LEHD-FMK conjugated to sulfo-rhodamine (Red-LEHD-FMK) as the fluorescent *in situ* 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 fluorescence plate reader.

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-LEHD-FMK | 100 μ L |
| Wash Buffer | 2 x 100 mL |
| Z-VAD-FMK | 100 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. Staining Procedure

- a) Induce apoptosis in cells (1×10^6 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 μ 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 μ l each of the induced and control cultures into eppendorf tubes.
- c) Add 1 μ l of Red-LEHD-FMK into each tube and incubate for 0.5-1 hour at 37°C incubator with 5% CO₂.
- 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 of the un-induced control.

2. Quantification by Flow Cytometry:

For flow cytometric analysis, re-suspend cells in 300 μ l of 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 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 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 μ l 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 Initiator Caspases

| Caspase | Cleavage motif | Inhibitor motif | Cross-reactivity with other caspase: | | | | | | | | | | | |
|-------------------|----------------|-----------------|--------------------------------------|---|---|---|---|---|---|---|----|----|--|---|
| | | | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | | |
| Caspase 2 | VDVAD | | | | Y | | | | | Y | | | | |
| Caspase 8 | IETD | IETD, LETD | | | Y | | | | Y | | | | | Y |
| Caspase 9 | LEHD | | | | Y | | | | Y | | Y | | | Y |
| Caspase 10 | AEVD | | | | Y | | | | | Y | Y? | | | |

6. 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 |

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “*contact us*” on www.abcam.com for the phone number for your region).

UK, EU and ROW

Email: technical@abcam.com | Tel: +44-(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com |
Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel:
01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com
Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 400 921 0189 / +86 21 2070 0500

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp