

ab65659

Caspase 4 Assay Kit (Colorimetric)

Instructions for Use

For the rapid, sensitive and accurate measurement of Caspase 4 activity in cell and tissue lysates

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

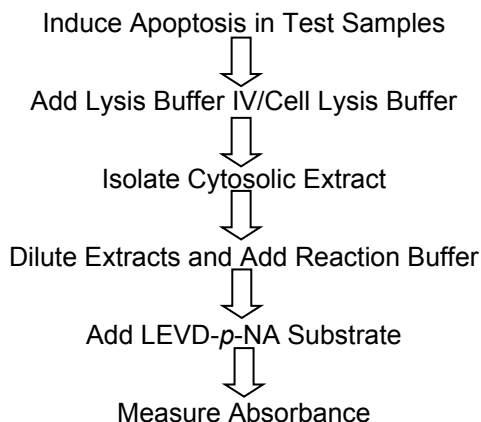
| | |
|--|---|
| Table of Contents | 1 |
| 1. Overview | 2 |
| 2. Protocol Summary | 2 |
| 3. Components and Storage | 3 |
| 4. Assay Protocol | 5 |
| 5. Data Analysis | 6 |
| 6. Factors to consider for caspase activity assays | 7 |
| 7. Troubleshooting | 9 |

1. Overview

Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells. Abcam's Caspase 4 Assay Kit (Colorimetric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence LEVD.

The assay is based on spectrophotometric detection of the chromophore *p*-nitroanilide (*p*-NA) after cleavage from the labeled substrate LEVD-*p*-NA. The *p*-NA light emission can be quantified using a spectrophotometer or a microtiter plate reader at 400 or 405nm. Comparison of the absorbance of *p*-NA from a treated sample with an untreated control allows determination of the fold increase in Caspase 4 activity.

2. Protocol Summary



3. Components and Storage

A. Kit Components

| Item | Quantity |
|---|-------------|
| Lysis Buffer IV/Cell Lysis Buffer | 100 mL |
| 2X Reaction Buffer I/2X Reaction Buffer | 4 x 2 mL |
| LEVD-pNA/LEVD-p-NA(4mM) | 500 μ L |
| DTT I/DTT(1M) | 400 μ L |
| Dilution Buffer II/Dilution Buffer | 100 mL |

* Store kit at -20°C. All reagents are stable for 1 year under proper storage conditions.

- Aliquot enough 2X Reaction Buffer I/2X Reaction Buffer for the number of assays to be performed. Add DTT I/DTT to the 2X Reaction Buffer I/2X Reaction Buffer immediately before use (10 mM final concentration: add 10 μ l of 1.0 M DTT I/DTT stock per 1ml of 2X Reaction Buffer I/2X Reaction Buffer).
- Protect LEVD-p-NA from light.

- After thawing, store Lysis Buffer IV/Cell Lysis Buffer, 2X Reaction Buffer I/2X Reaction Buffer, and Dilution Buffer II/Dilution Buffer at +4°C.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Colorimetric microplate reader or spectrophotometer
- 96-well plate
- Micro-quartz and regular cuvettes
- Orbital shaker

4. Assay Protocol

1. Treat samples with the desired method to induce caspase activity. TNF alpha is a common treatment to induce inflammation (and Caspase 4). Concurrently, incubate a separate culture without treatment to use as a negative control.

Note:

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

2. Count cells and pellet 5×10^6 cells.
3. Re-suspend in 50 μ l of chilled Lysis Buffer IV/Cell Lysis Buffer and incubate on ice for 10 min.
4. Centrifuge for 1min in a microcentrifuge (10,000 x g).
5. Transfer supernatant (cytosolic extract) to a fresh tube and keep on ice.
6. Assay protein concentration.
7. Dilute 200-300 μ g protein to 50 μ l Lysis Buffer IV/Cell Lysis Buffer for each assay.
8. Add 50 μ l of 2X Reaction Buffer I/2X Reaction Buffer (containing 10 mM DTT I/DTT) to each sample. Add 5 μ l of the LEVD-

pNDA/4 mM LEVD-*p*-NA substrate (200 μ M final conc.). Incubate at 37°C for 1-2 hours.

9. Read samples at 400 or 405 nm in a microtiter plate reader, or spectrophotometer using a 100 μ l micro quartz cuvette, or dilute sample to 1 ml with Dilution Buffer II/Dilution Buffer and using regular cuvette (Dilution of the samples proportionally decreases the reading).

You may also perform the entire assay in a 96-well plate.

5. Data Analysis

Fold-increase in LEVD-dependent Caspase activity can be determined by comparing the results of induced samples with the level of the un-induced control.

Note:

Background reading from cell lysates and buffers should be subtracted from the readings of both induced and the un-induced samples before calculating fold increase in Caspase 4 activity.

6. 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, 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.

Inflammatory Caspases

| Caspase | Cleavage motif | Inhibitor motif | Cross-reactivity with other caspase: | | | | | | | | | | | |
|------------|----------------|-----------------|--------------------------------------|---|---|--------|---|---|---|---|---|----|--|--|
| | | | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | | |
| Caspase 1 | YVAD | | | | | Y ? | Y | | | | | | | |
| Caspase 4 | LEVD | LEHD* | | | | | Y | | | | | | | |
| Caspase 5 | WEHD | LEHD* | Y | | | Y | | | | | | | | |
| Caspase 12 | ATAD | | | | | | | | | | | | | |

** inhibits at high concentration*

7. Troubleshooting

| Problem | Reason | Solution |
|-------------------|--|---|
| Assay not working | Cells did not lyse completely | Re-suspend the cell pellet in the lysis buffer and incubate as described in the datasheet |
| | Experiment was not performed at optimal time after apoptosis induction | Perform a time-course induction experiment for apoptosis |
| | Plate read at incorrect wavelength | Ensure you are using appropriate reader and filter settings (refer to datasheet) |
| | Old DTT used | Always use freshly thawed DTT in the cell lysis buffer |
| High Background | Increased amount of cell lysate used | Refer to datasheet and use the suggested cell number to prepare lysates |
| | Increased amounts of components added due to incorrect pipetting | Use calibrated pipettes |
| | Incubation of cell samples for extended periods | Refer to datasheet and incubate for exact times |
| | Use of expired kit or improperly stored reagents | Always check the expiry date and store the individual components appropriately |
| | Contaminated cells | Check for bacteria/ yeast/ mycoplasma contamination |

| Problem | Reason | Solution |
|-------------------------------|---|---|
| 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 datasheet for appropriate cell number |
| | Use of samples stored for a long time | Use fresh samples or aliquot and store and use within one month for the assay |
| | Incorrect setting of the equipment used to read samples | Refer to datasheet and use the recommended filter setting |
| | Allowing the reagents to sit for extended times on ice | Always thaw and prepare fresh reaction mix before use |
| Samples with erratic readings | Uneven number of cells seeded in the wells | Seed only equal number of healthy cells (correct passage number) |
| | Samples prepared in a different buffer | Use the cell lysis buffer provided in the kit |
| | Adherent cells dislodged and lost at the time of experiment | Perform experiment gently and in duplicates/triplicates; apoptotic cells may become floaters |
| | Cell/ tissue samples were not completely homogenized | Use Dounce homogenizer (increase the number of strokes); observe efficiency of lysis under microscope |
| | Samples used after multiple freeze-thaw cycles | Aliquot and freeze samples, if needed to use multiple times |
| | Presence of interfering substance in the sample | Troubleshoot as needed |
| | Use of old or inappropriately stored samples | Use fresh samples or store at correct temperatures until use |

| | | |
|--------------------|---|---|
| Unexpected results | Measured at incorrect wavelength | Check the equipment and the filter setting |
| | Cell samples contain interfering substances | Troubleshoot if it interferes with the kit (run proper controls) |
| General Issues | Improperly thawed components | Thaw all components completely and mix gently before use |
| | Incorrect incubation times or temperatures | Refer to datasheet & verify the correct incubation times and temperatures |
| | Incorrect volumes used | Use calibrated pipettes and aliquot correctly |
| | Air bubbles formed in the well/tube | Pipette gently against the wall of the well/tubes |
| | Substituting reagents from older kits/ lots | Use fresh components from the same kit |
| | Use of a different 96-well plate | Fluorescence: Black plates; Absorbance: Clear plates |

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).

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

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)