



**Caspase-9/Mch6/Apaf-3
Colorimetric Protease Assay
Catalog# KHZ0101 (25 Tests)
KHZ0102 (200 Tests)**



www.invitrogen.com
Invitrogen Corporation
542 Flynn Rd, Camarillo, CA 93012
Tel: 800-955-6288
E-mail: techsupport@invitrogen.com

INTENDED USE

The *ApoTarget*[™] Caspase-9 Protease Assay is to be used for the *in vitro* determination of Caspase-9 proteolytic activity in lysates of mammalian cells. This test is designed for research use only. It is not to be used in diagnostic procedures.

PRINCIPLE OF THE METHOD

Caspase-9 (also known as Mch6, Apaf-3 or ICE-LAP-6) is a member of the Interleukin-1 β Converting Enzyme (ICE) family of cysteine proteases. Caspase-9 exists in cells as an inactive 46 kDa proenzyme, called pro-Caspase-9. Pro-Caspase-9 is cleaved to active 35 and 10 kDa subunits by complexing with a cofactor, Apaf-1, through a caspase recruitment domain (CARD) and with cytochrome c during apoptosis. The downstream substrates of Caspase-9 include Caspases-3 and -7. Caspase-9 plays a critical role in the survival of neuronal progenitors within the embryonic forebrain and cortex. Phosphorylation of Caspase-9 by Akt, a serine/threonine kinase, inhibits its protease activity.

The Caspase-9 colorimetric protease assay provides a simple and convenient means for quantitating the enzyme activity of caspases that recognize the amino acid sequence, LEHD. The kit includes substrate and optimized buffers. The substrate, LEHD-*p*NA, is composed of the chromophore, *p*-nitroanilide (*p*NA), and a synthetic tetrapeptide, LEHD (Leu-Glu-His-Asp). Upon cleavage of the substrate by Caspase-9 or related Caspases, absorption of light by free *p*NA can be quantified using a spectrophotometer or a microtiter plate reader at 400 or 405 nm. Comparison of the absorbance of *p*NA from apoptotic sample with an uninduced control allows determination of the fold increase in Caspase-9 activity.

REAGENTS PROVIDED

Note: Store kit at -20°C. Once opened, store Cell Lysis Buffer, 2x Reaction Buffer, and Dilution Buffer at 2-8°C. Refer to reagent label for expiration.

1. Cell Lysis Buffer (25 mL or 100 mL): Tris buffered saline containing detergent.
2. 2x Reaction Buffer (2 mL or 16 mL): contains buffered saline, glycerol and detergent.
3. Substrate (125 μ L or 1000 μ L): contains 4 mM of synthetic peptide (sequence: LEHD) conjugated to the chromophore, *p*NA, in DMSO.
4. Dilution Buffer (25 mL or 200 mL): buffered solution containing HEPES, Tris and NaCl.
5. DTT (100 μ L or 800 μ L): contains 1 M dithiothreitol.

MATERIALS NOT PROVIDED

1. Spectrophotometer (and cuvettes) or microplate reader capable of measurement at 400-405 nm.
2. Calibrated adjustable precision pipettes, preferably with disposable plastic tips.
3. Protein measurement method such as Bradford protein assay.
4. Tubes appropriate for holding cells during induction of apoptosis.
5. Microcentrifuge.
6. Reaction tubes or 96-well microplate.

PROCEDURAL NOTES/LAB QUALITY CONTROL

1. When not in use, kit components should be stored refrigerated or frozen as indicated on vial or bottle labels. All reagents should be warmed to room temperature before use.
2. If not analyzed immediately, samples should be stored at -20°C or lower.
3. It is recommended that all samples and controls be run in duplicate.
4. Protect LEHD-*p*NA from light.
5. Cover or cap all reagents when not in use.
6. Do not mix or interchange different reagent lots from various kits.

7. Do not use reagents beyond the expiration date of the kit.
8. Set negative controls using same number of cells or same amount of lysates without adding LEHD-*p*NA substrate.

WARNINGS AND PRECAUTIONS

1. This kit is intended for research use only. It is not to be used for diagnostic procedures.
2. Never pipette by mouth.
3. Do not eat, drink or smoke in the laboratory areas. All blood components and biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

ASSAY PROCEDURE

1. Induce apoptosis in cells by desired method. Concurrently incubate a control culture without induction.
2. Count cells and pellet $3\text{--}5 \times 10^6$ cells per sample.
3. Resuspend cells in 50 μL of chilled Cell Lysis Buffer and incubate cells on ice for 10 minutes.
4. Centrifuge for 1 minute in a microcentrifuge (10,000 \times g).

Note: This step is not recommended if using the fluorometric caspase assay.

5. Transfer supernatant (cytosol extract) to a fresh tube and put on ice.
6. Assay protein concentration by any standard method.
7. Dilute each cytosol extract to a concentration of 50-200 μg protein per 50 μL Cell Lysis Buffer (1-4 mg/mL).
8. Determine the number of samples to be measured and aliquot enough 2x Reaction Buffer into a glass tube (assuming 50 μL of 2x Reaction Buffer per sample). Add DTT to the 2x Reaction Buffer immediately before use (10 mM final concentration: add 10 μL of 1.0 M DTT stock per 1 mL of 2x Reaction Buffer).
9. Add 50 μL of 2x Reaction Buffer (containing 10 mM DTT) to each sample.
10. Add 5 μL of the 4 mM LEHD-*p*NA substrate (200 μM final concentration) and incubate at 37°C for 2 hours. Keep the samples in the dark during incubation.
11. Read samples at 400 nm or 405 nm in a microplate reader, or spectrophotometer using a 100 μL micro-quartz cuvette, or dilute sample to 1 mL with Dilution Buffer and use a regular cuvette.

Note: Dilution of the samples proportionally decreases the optical density. You may also perform the entire assay directly in a 96-well plate.

12. Fold-increase in Caspase-9 activity should be determined by direct comparison to the level of the uninduced control.

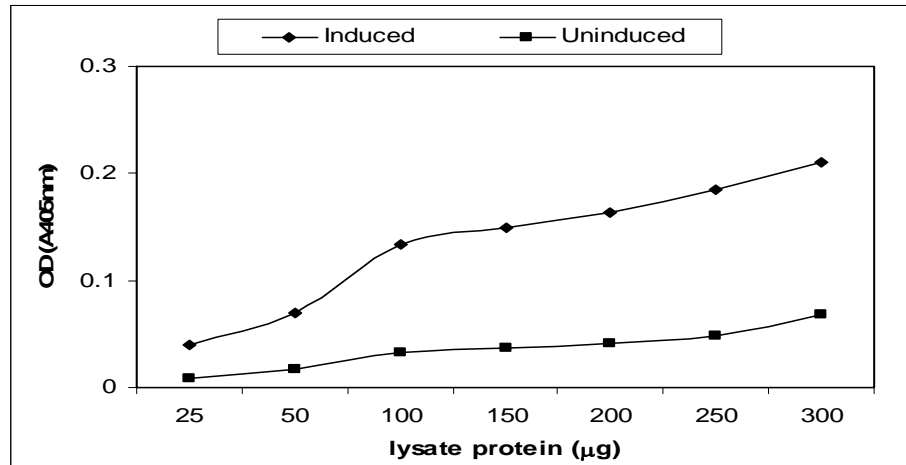
Note: Background absorbances from cell lysates and buffers should be subtracted from the absorbances of both induced and the uninduced samples before calculating fold-increase in Caspase-9 activity.

TYPICAL DATA

The following absorbance data were obtained from Jurkat cells. Apoptosis was induced by incubating Jurkat cells with 10 $\mu\text{g/mL}$ Camptothecin for 6 hours. The assay was performed according to the procedure described above.

LIMITATIONS OF THE PROCEDURE

This kit provides a simple and convenient method to detect Caspase-9 activity of apoptotic cells. A relatively high concentration of DTT (10 mM) is required for full activity of Caspase enzymes. Make sure that DTT is added to Reaction Buffer when the assay is carried out. Otherwise, unexpected low Caspase-9 activity will occur. Turbidity, lipids or particulate material in samples can decrease the assay precision.



REFERENCES:

1. Zou, H., et al. (1999) An APAF-1 cytochrome c multimeric complex is a functional apoptosome that activates procaspase-9. *J. Biol. Chem.* 274:11549-11556.
2. Slee, E.A., et al. (1999) Ordering the cytochrome c-initiated caspase cascade: hierarchical activation of caspases-2, -3, -6, -7, -8, and -10 in a caspase-9-dependent manner. *J. Cell. Biol.* 144(2):281-292.
3. Hakem, R., et al. (1998) Differential requirement for caspase-9 in apoptotic pathways in vivo. *Cell* 94(3):339-352.
4. Kuida, K., et al. (1998) Reduced apoptosis and cytochrome c-mediated caspase activation in mice lacking caspase-9. *Cell* 94(3):325-337.
5. Chou, J.J., et al. (1998) Solution structure of the RAIDD CARD and model for CARD/CARD interaction in caspase-2 and caspase-9 recruitment. *Cell* 94(2):171-180.
6. Thornberry, N.A., et al. (1997) A combinatorial approach defines specificities of members of the caspase family and granzyme B. Functional relationships established for key mediators of apoptosis. *J. Biol. Chem.* 272(29):17907-17911.
7. Talanian, R.V., et al. (1997) Substrate specificities of caspase family proteases. *J. Biol. Chem.* 272(15):9677-9682.
8. Stennicke, H.R., et al. (1997) Biochemical characteristics of Caspases-3, -6, -7, and -8. *J. Biol. Chem.* 272(41):25719-25723.

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.