



CV-Caspase 3&7 detection kit

Catalog # BML-AK118

96 Well Plate Kit



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON- INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

TABLE OF CONTENTS



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

Introduction.....	4
Materials Provided.....	6
Storage.....	6
Other Materials Needed.....	6
Detection Equipment	7
Protocol	8
References	22
Contact Information	24

INTRODUCTION

Enzo Life Sciences' CV-Caspase 3&7 assay kit enables researchers to detect and monitor apoptosis over time via intracellular caspase activity *in vitro*. The CR(DEVD)₂ reagent is a non-cytotoxic substrate that fluoresces upon cleavage by active caspase 3&7 enzymes. It measures the intracellular process of apoptosis instead of a side-effect, such as the turnover of phosphatidyl serine.

To use CR(DEVD)₂, add the reagent directly to the cell culture media, incubate, and analyze. Because CR(DEVD)₂ is cell-permeant, it easily penetrates the cell membrane and the membranes of the internal cellular organelles - no lysis or permeabilization steps are required. CR(DEVD)₂ will enter the cell in a non-fluorescent state. If caspase 3&7 enzymes (DEVDases) are active, they will cleave CR(DEVD)₂ and release the cresyl violet fluorophore, which will begin to fluoresce red. The red fluorescent product will stay inside the cell and will often aggregate inside lysosomes (Figures 1 and 3); caspases are not lysosomal. As protease activity progresses and more CR(DEVD)₂ substrate is cleaved, the red fluorescent signal will intensify, enabling researchers to watch the color develop over time (Figure 2) and quantify apoptosis (Figures 6 and 9).

There is no interference from pro-caspases nor inactive forms of the enzymes. If the treatment is causing cell death via apoptosis, apoptotic cells will have elevated levels of caspase 3&7 activity relative to non-apoptotic or negative control cells.

Up-regulation and initiation of the caspase enzyme cascade is the central driving force behind apoptosis³. Although a number of other intracellular enzyme families, including the cathepsins, calpains, and granzymes, participate in the cell break-down mechanism, the caspase cascade occupies the central effector role in the cell suicide process^{21,25,8,12}. Like other intracellular proteases, caspases are initially synthesized as inactive zymogen precursors that can be rapidly activated upon auto and heterologous enzymatic processing at specific sites containing an aspartic acid¹⁷. Caspase 3 is the predominant effector caspase in apoptosis (with few exceptions, such as MCF-7 cells, which are deficient in caspase 3; Figures 1, 3, and 4).

Caspase enzymes cleave proteins and are classified as cysteine proteases based on the mechanism of substrate hydrolysis at their active site. Caspases specifically recognize a 3 or 4 amino acid sequence which must include an aspartic acid residue (D)

in the P1 position. This C-terminal residue is the target for the cleavage reaction at the carbonyl end¹⁵. Enzo Life Sciences' CV-caspase 3&7 probe contains a 4 amino acid sequence, aspartylglutamylvalanyl aspartic acid (DEVD), which is the optimal target sequence for caspase 3 as well as caspase 7²². Two copies of this sequence are coupled to a photostable red fluorophore, cresyl violet, to create the CV-caspase 3&7 substrate CR(DEVD)₂. The intact CR(DEVD)₂ substrate reagent is not fluorescent; the fluorescent signal is generated by the cresyl violet fluorophore once CR(DEVD)₂ is cleaved by caspase 3 or caspase 7. CR(DEVD)₂ has been shown to work in human, rat, and mouse cells, among other species. A baseline level of DEVDase activity is present in all cell lines. Apoptotic cells will fluoresce red and have pronounced red lysosomes and mitochondria. Healthy and nonapoptotic cells will exhibit very low levels of background red fluorescence evenly distributed throughout the cell (Figures 1-3, 6-8). This background level of substrate activity could be the result of constitutively synthesized serine proteases that target analogous aspartic acid sequences for hydrolysis. Cells in more advanced stages of apoptosis, containing peak levels of DEVDase activity, will appear brighter red than cells in earlier stages of apoptosis.

The fluorophore, cresyl violet, fluoresces red when excited at 550-590 nm²³. The red fluorescent signal can be monitored with a fluorescence microscope (Figures 1-3, 6-8) or plate reader (Figure 9). It has an optimal excitation of 592nm and emission of 628 nm². At these higher excitation wavelengths, the amount of cell-mediated auto-fluorescence is minimal²². The excitation peak is rather broad allowing good excitation efficiency at 540-560 nm. Hoechst stain is included with the kit to concurrently label nuclei after labeling with CR(DEVD)₂ (Figures 1 and 6). Acridine orange (AO) is also included in the kit to identify lysosomes and other intracellular organelles (Figures 4 and 5).

MATERIALS PROVIDED

1. 1 vial of CV-Caspase 3&7 Substrate Reagent (CR(DEVD)₂), approximately 25 tests, BML-KI224
2. 1 vial of Hoechst 33342, 200 µg/mL (1 mL), BML-KI189
3. 1 vial of Acridine Orange, 1 mM (0.5 mL), BML-KI223

STORAGE

Store the unopened kit and each unopened component at 2-8°C until the expiration date. Once reconstituted with DMSO, use CR(DEVD)₂ immediately, or store at ≤-20°C up to 6 months, protected from light and thawed no more than twice during that time.

OTHER MATERIALS NEEDED

- DMSO, 100-400 µL to reconstitute CR(DEVD)₂
- diH₂O, 400-1600 µL to dilute CR(DEVD)₂
- Phosphate buffered saline (PBS) pH 7.4, 100 mL
- Cultured cells treated with the experimental conditions ready for staining
- Reagents to induce apoptosis and create controls, such as staurosporine (BML-EI156) or camptothecin (ALX-350-015)
- Hemocytometer
- Centrifuge at <200g
- 15 mL polystyrene centrifuge tubes
- Sterile black 96-well microtiter tissue culture plates, round or flat bottom.
- Slides and coverslips
- Ice or refrigerator

DETECTION EQUIPMENT

CR(DEVD)₂ excites at 550 nm and emits >610 nm. It has an optimal excitation and emission wavelength tandem of 592 nm and 628 nm, respectively. Hoechst stain can be visualized using a UV-filter with excitation at 365 nm and emission at 480 nm (Section 5). AO excites at 480 nm and emits >540 nm (Section 6). Select a filter combination that best approximates these settings.

Fluorescence microscope: Use an excitation filter of 550 nm (540-560 nm) and a long pass >610 nm emission/barrier filter pairing. Using this excitation/emission filter pairing, cells should stain red with brightly stained vacuoles and lysosomes (Figures 2 and 3). If the samples were stained with both CR(DEVD)₂ and Hoechst, the dual staining properties can be examined using a multi-wavelength filter (Figures 1 and 6). Hoechst stain can be seen using a UV-filter with excitation at 365 nm and emission at 480 nm. As AO exhibits a very broad emission range, one of several filter pairings may be used. The same excitation/emission filter pairings used to view CR(DEVD)₂ may be used for AO: a 550 nm (540 – 560 nm) excitation filter with a long pass >610 nm emission/barrier filter. With this pairing, lysosomes appear red. When illuminating AO with a blue light (480 nm) excitation filter, a green light (540-550 nm) emission/barrier filter combination works well. Lysosomes will appear yellowish green instead of red (Figures 4 and 5).

Fluorescence plate reader: CR(DEVD)₂ has an optimal excitation and emission wavelength tandem of 592 nm and 628 nm, respectively. Use a fluorescence plate reader with excitation at 590 nm and emission at 630-640 nm. If available, use a cut-off filter at 630 nm to filter out shorter wavelength excitation interference (Figure 9).

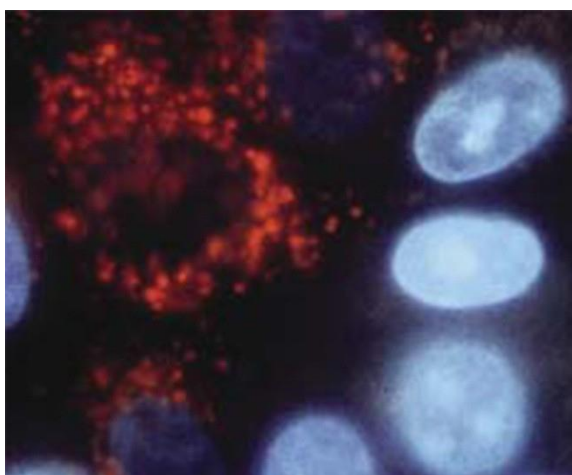
PROTOCOL

1. Experimental Preparation

Staining apoptotic cells with CR(DEVD)₂ can be completed within a few hours. However, CR(DEVD)₂ is used with living cells, which require periodic maintenance and cultivation several days in advance. In addition, once the proper number of cells has been cultivated, time must be allotted for the experimental procedure or apoptosis induction (which typically requires a 3-6 hour incubation at 37°C).

As CR(DEVD)₂ detects caspase-mediated apoptosis, plan the experiment so that it will be diluted and administered at the time when caspases are expected to be activated in the cells. The recommended volume of the CR(DEVD)₂ staining solution is 0-20 µL per 300 µL of cells at 10⁶ cells/mL, but the ideal amount may vary based on the experimental conditions and method of analysis. Culture cells to a density optimal for the specific experiment or apoptosis induction protocol. Cell density should not exceed 10⁶ cells/mL as cells cultivated in excess of this concentration may begin to naturally enter apoptosis due to nutrient deprivation or the accumulation of cell degradation products in the media.

FIGURE 1: DUAL STAINING OF MCF-7 CELLS



Apoptotic MCF-7 cells were dually stained using Enzo Life Sciences' CR(DEVD)₂ fluorogenic caspase 3&7 substrate and Hoechst 33342 nuclear stain (BML-AK118-0001). MCF-7 cells were exposed to 0.15 µM camptothecin (ALX-350-015) for 24 hours at 37°C, then stained with CR(DEVD)₂ for 30 minutes at 37°C, washed twice in PBS, and supravivally stained with 1 µg/mL Hoechst stain for about 10 minutes. Using the Nikon Microphot FXA system with multi-wavelength filter pairs (UV for Hoechst stain and green light for CR(DEVD)₂), apoptotic cells bearing orange-red lysosomal bodies with less intense blue nuclei are intermixed with non-apoptotic cells bearing bright blue nuclei and absent or reduced orange-red lysosomal staining. Photo provided by Dr. Zbigniew Darzynkiewicz at Brander Cancer Research Center Institute, New York City, NY.

Cells undergoing apoptosis will generate a stronger red fluorescence with CR(DEVD)₂ than non-apoptotic cells of the same lineage. To optimize this assay, determine the greatest difference in the fluorescent signal between induced and non-induced cell populations. Adjust the amount of CR(DEVD)₂ substrate used to stain cells and the incubation time. Hoechst stain can be used with CR(DEVD)₂ to label nuclei (Figures 1 and 6). Because of the overlap in emissions, dual staining of cells with both CR(DEVD)₂ and AO will yield confusing results and is not recommended; these dyes should be used separately. Do not use CR(DEVD)₂ with paraffin-embedded tissues as the chemicals used for paraffin-embedding may denature and inactivate the substrate.

2. Controls

It is highly recommended that two sets of controls be run: one positive population of cells that was induced to undergo apoptosis or trigger caspase 3/7 activity; and a placebo population of cells that received just the vehicle used to deliver the apoptosis-inducing agent (Figure 6). Create positive controls by inducing apoptosis (Section 3). Create negative controls by culturing an equal volume of non-induced cells for every labeling condition. The negative control and induced positive control populations should contain similar quantities of cells. For example, if labeling with CR(DEVD)₂, Hoechst stain, and Acridine Orange, make 10 control populations:

- a. Unlabeled, induced and non-induced populations.
- b. CR(DEVD)₂-labeled, induced and non-induced populations.
- c. CR(DEVD)₂- and Hoechst-labeled, induced and non-induced populations.
- d. Hoechst-labeled, induced and non-induced populations.
- e. AO-labeled, induced and non-induced populations.

FIGURE 2: WATCH CASPASE ACTIVITY IN REAL TIME



REC:MYC immortalized rat fibroblasts were seeded in a 12-well plate at 1×10^4 cells/mL and irradiated (4 Gy, X-ray) the following day. Enzo Life Sciences' Caspase 3&7 CR(DEVD)₂ Substrate Reagent (BML-KI224) was added, and cells were photographed for 16 hours using an inverted Nikon TE2000 microscope with a CCD camera from Hamamatsu and PCI software from Compix. The red fluorescence became brighter as caspase activity and apoptosis progressed. Data courtesy of Dr. Martin Purschke, Massachusetts General Hospital.

3. Apoptosis Induction

CR(DEVD)₂ works with your current apoptosis protocols - induce apoptosis or trigger DEVDase activity as you normally would, then label cells with CR(DEVD)₂. For example, apoptosis may be induced with:

- a. 2-4 μ g/mL camptothecin for >4 hours
- b. 1-2 μ M staurosporine for >4 hours

4. CR(DEVD)₂ Substrate Reagent

CR(DEVD)₂ is supplied as a lyophilized powder that may be slightly visible as an iridescent sheen inside the vial. It must first be reconstituted in DMSO, forming the stock concentrate, and then diluted 1:5 in diH₂O to form the final staining solution. For best results, the staining solution should be prepared immediately prior to use. However, the stock concentrate may be stored at $\leq 20^\circ\text{C}$ for future use. Protect from light and use gloves when handling.

1. Create the stock solution by reconstituting CR(DEVD)₂.
 - a) Use 100 μ L DMSO to reconstitute the vial BML-KI224.
2. Gently vortex or swirl the vial, allowing the DMSO to travel around the base of the vial until completely dissolved. At room temperature (RT), this should take just a few minutes. The stock solution should appear red. Once reconstituted, it may be stored at $\leq -20^\circ\text{C}$ up to 6 months protected from light and thawed no more than twice during that time. If using immediately, dilute

in diH₂O to form the staining solution. If not diluting within 1 hour, aliquot and freeze.

3. Immediately prior to staining the samples, dilute the stock solution 1:5 in diH₂O to form the staining solution. Use the staining solution within 15 minutes of dilution to prevent substrate hydrolysis.
 - b) Add 400 µL diH₂O to dilute the vial BML-KI224. The vial contains 100 µL of the stock (1a); this yields 500 µL of the staining solution. Use immediately.
 - c) For other amounts, dilute the stock 1:5 in diH₂O. For example, add 10 µL stock to 40 µL diH₂O; this yields 50 µL of the staining solution. Use immediately.
 - d) Mix by inverting or vortexing the vial at RT.

5. Hoechst Stain

Hoechst 33342 is a cell-permeant nuclear stain that emits blue fluorescence when bound to double stranded DNA. It is used: a) to stain the nuclei of living or fixed cells; b) to distinguish condensed pyknotic nuclei in apoptotic cells; and c) for cell cycle studies. When bound to nucleic acids, the maximum absorption is 350 nm and the maximum emission is 480 nm. It is revealed under a microscope using a UV-filter with excitation at 365 nm and emission at 480 nm. Hoechst Stain (BML-KI189) is provided ready- to-use at 200 µg/mL. Hoechst Stain can be used concurrently with the CR(DEVD)₂ to label nuclei (Figures 1 and 6).

Warning: Hoechst Stain is a potential mutagen. Gloves, protective clothing, and eye wear are strongly recommended. When disposing, flush sink with copious amounts of water. See MSDS for further information.

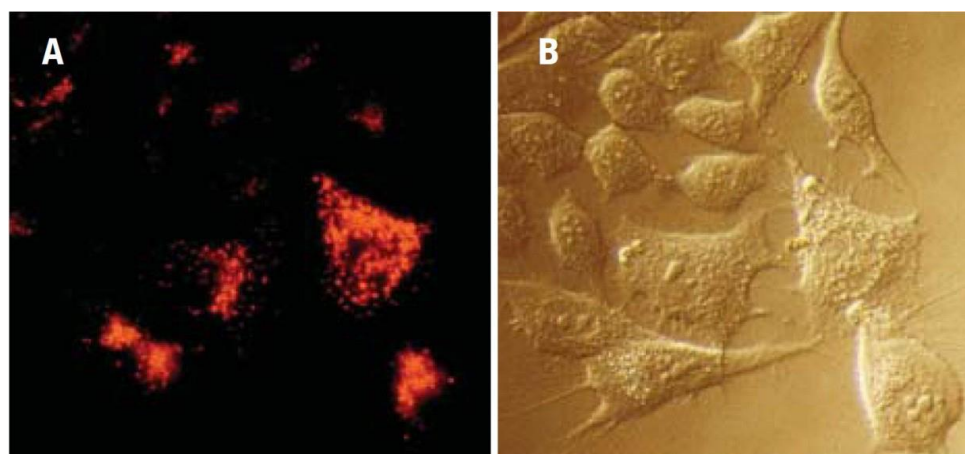
6. Acridine Orange

Acridine Orange (AO) is a chelating dye and can be used to reveal lysosomes, nuclei, and nucleoli (Figures 4 and 5). 0.5 mL of AO is provided at 1 mM (ML-KI223). AO may be used neat or diluted in diH₂O or media prior to pipetting into the cell suspension. Always protect AO from bright light.

Lysosomal structures can be visualized by staining with AO at 0.5- 5.0 µM. This concentration range can be obtained by diluting the AO reagent stock 1:2,000-1:200 (0.05-0.5% v/v)

into the final cell suspension. For example, if using AO at 1.0 μM in the final cell suspension, it must be diluted 1:1,000. First dilute it 1:100 in PBS; e.g., put 10 μL AO into 990 μL PBS. Pipette the diluted AO into the cell suspension at approximately 1:10; e.g., put 50 μL diluted AO into 450 μL cell suspension.

FIGURE 3: CASPASE ACTIVITY IN MCF-7 CELLS



DEVDase activity was detected in apoptotic MCF-7 cells using Enzo Life Sciences' CR(DEVD)₂ caspase 3&7 fluorogenic substrate. Apoptosis was induced in MCF-7 cells by treating them with 0.15 μM camptothecin (ALX-350-015) for 24 hours at 37°C. Cells were then exposed to CR(DEVD)₂ for 60 minutes at 37°C. DEVDase activity is demonstrated by the appearance of orange-red lysosomal bodies within the cytoplasm of the cell (A). The photograph was taken on a Nikon Microphot FXA system at 541-551 nm excitation with a long pass >640 nm barrier filter (compare with AO staining in Figure 4). Photo B shows the corresponding interference contrast image of the cells. Data courtesy of Dr. Zbigniew Darzynkiewicz (Brander Cancer Research Center Institute, New York City, NY).

As AO exhibits a very broad emission range, one of several filter pairings on the fluorescence microscope can be used to view this stain. The same excitation/emission filter pairings used to view CR(DEVD)₂ may be used: an excitation filter of 550 nm (540-560 nm) and a long pass >610 nm emission/barrier filter pair. With this pairing, the lysosomes appear red instead of yellowish green.

When illuminating with a blue light (480 nm) excitation filter, a green light (540-550 nm) emission/barrier filter combination works well. Lysosomes will appear yellowish green. As this filter combination is very close to the maximum emission of AO, the slide may appear too bright. Excess AO may be removed by washing cells prior to viewing.

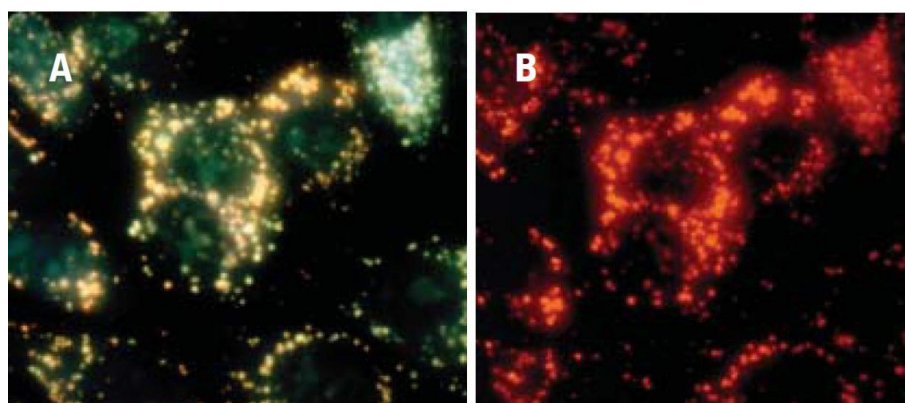
Because of the overlap in emissions, dual staining of cells with both CR(DEVD)₂ and AO will yield confusing results. Therefore, these dyes should be used separately. Warning:

AO is a potent mutagen and probable carcinogen. Use gloves, protective clothing, and eye wear. When disposing, flush sink with copious amounts of water. See MSDS for further information.

7. Microscopy Analysis of Suspension Cells

1. Prepare cell populations. Initial cell concentrations should be $3-5 \times 10^5$ cells/mL and should not exceed 7×10^5 cells/mL, as cells cultivated in excess of this concentration may begin to naturally enter apoptosis.
2. Expose cells to the experimental conditions and create positive and negative controls (Section 2) or induce cells to undergo apoptosis (Section 3).
3. When ready to label with the staining solution, cell concentrations should be $1-2 \times 10^6$ cells/mL for best viewing. Density can be determined by counting cell populations on a hemocytometer. If necessary, concentrate cells by gentle centrifugation at $200 \times g$ for 3-8 minutes. Remove the supernatant and resuspend with cell culture media or PBS.
4. Transfer 300 μ L into 12 x 75 mm glass or polypropylene tubes. If desired, larger cell volumes can be used, but additional CR(DEVD)₂ staining solution may be required.

FIGURE 4: ACRIDINE ORANGE STAINING OF MCF-7 CELLS

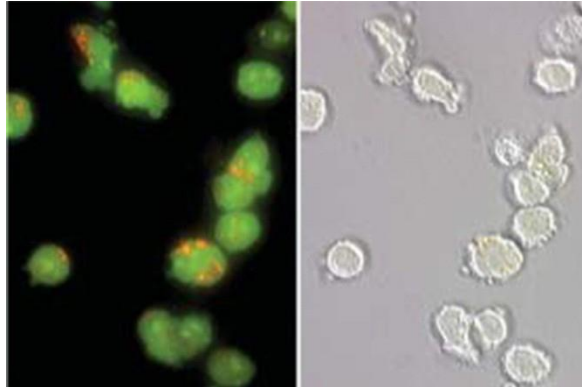


Apoptosis was induced in MCF-7 cells by treating them with $0.15 \mu\text{M}$ camptothecin (ALX-350-015) for 24 hours at 37°C . Cells were stained with AO in PBS for 30 minutes, then washed twice in PBS. Cells were photographed with a Nikon Microphot-FXA epifluorescence microscope at 40X using either a blue light excitation (492 nm) with a 540-550 nm emission filter (A, lysosomes appear yellowish green), or green light excitation (540 nm) with a long pass >640 nm barrier filter (B, lysosomes appear red; compare with CR(DEVD)₂ staining of MCF cells in Figure 3, and

with Jurkat cells in Figure 5). Experiment performed in the laboratory of Dr. Zbigniew Darzynkiewicz (Brander Cancer Research Center Institute, New York City, NY).

5. Reconstitute CR(DEVD)₂ to form the stock solution (Section 4):
 - a. Use 100 μ L DMSO to reconstitute the vial (BML-KI224).
6. When ready to stain cells, dilute the stock 1:5 in diH₂O to form the CR(DEVD)₂ staining solution (Section 4):
 - a. Add 400 μ L diH₂O to dilute the vial (BML-KI224).
7. Add 10 μ L of the staining solution to each 300 μ L cell suspension and mix thoroughly. If different cell volumes are used, add the CR(DEVD)₂ staining solution at a ratio of approximately 1:30. For example, add 35 μ L CR(DEVD)₂ staining solution to 1,000 μ L of cell suspension forming a final volume of 1,035 μ L. Do not add CR(DEVD)₂ to cells that are to be labeled with AO; add a placebo instead, such as diH₂O (Step 10).
8. Incubate cells for 1 hour at 37°C and protect from light. Cells may settle on the bottom of the tubes; gently resuspend them by swirling cells every 20 minutes during the incubation to ensure even distribution of CR(DEVD)₂ substrate. After the incubation, cells can be stained with Hoechst Stain (Section 9), or unstained cells may be labeled with AO (Section 10).
9. If cells are to be labeled with Hoechst Stain, add it at approximately 0.5% v/v. Add 1.55 μ L Hoechst to 310 μ L cell suspension. Incubate 5 minutes at 37°C. Go to Step 11.
10. Because of the overlap in emissions, dual staining of cells with both CR(DEVD)₂ and AO is not recommended; the dyes should be used separately. To stain cells with AO:
 - a. Dilute AO to 1:2,000-1:200 (which is 0.05-0.5% v/v) into the final cell suspension. For example, if using AO at 1.0 μ M in the final cell suspension, first dilute it 1:100 in PBS; e.g., put 10 μ L AO into 990 μ L PBS. Pipette the diluted AO into the cell suspension at 1:10; e.g., add 35 μ L to 315 μ L cell suspension.
 - b. Incubate 30 minutes at 37°C.

FIGURE 5: ACRIDINE ORANGE STAINING OF JURKAT CELLS



Jurkat cells were stained with Acridine Orange (AO) in PBS for 60 minutes at 37°C. Jurkat cells stained with AO show orange lysosomal staining (photo A). Photomicrographs were taken using a Nikon Eclipse E800 photomicroscope using a 460- 500 nm excitation filter and a 505- 560 nm emission / barrier filter set at 300X. AO-stained lysosomes appear in photo A; photo B shows the corresponding DIC image of the cells (compare with Figure 4).

c. If viewing under the same filters used for CR(DEVD)₂ (excitation at 550 nm; emission >610 nm), cells may be viewed immediately after staining without a wash step - go to Step 11.

d. If viewing under blue (480 nm) excitation and green (540-550 nm) emission wavelengths, any excess AO may have to be washed away as the cells may appear too bright at this range. Brightness will depend on the type of microscope used and the cell line. To wash cells:

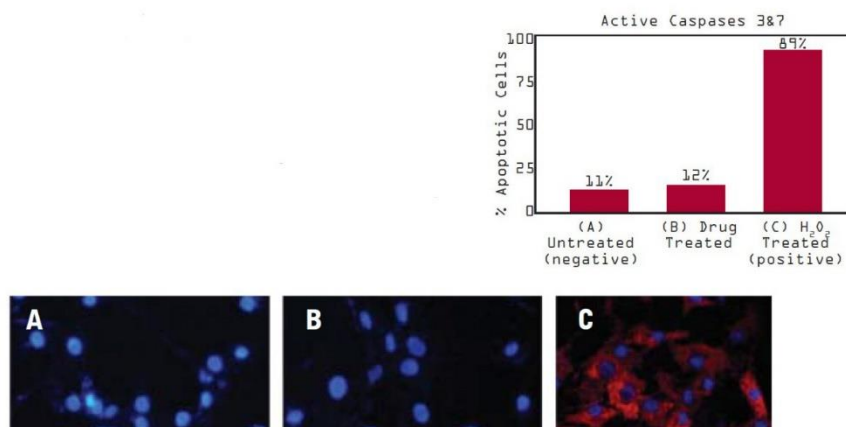
- i) Gently pellet cells at 200 x g for 3-8 minutes at RT.
 - ii) Remove and discard supernatant.
 - iii) Resuspend cells in 300 µL or a similar volume of PBS in which the cells were originally suspended.
11. Place 15-20 µL of cell suspension onto a microscope slide and cover with a coverslip.
 12. Observe cells using a fluorescence microscope equipped with an excitation filter of 550 nm (540- 560 nm) and a long pass >610 nm emission/barrier filter pairing. Select a filter combination that best approximates these settings. Using these filters, positive cells will appear red with brightly stained vacuoles and lysosomes.

If the samples were stained with both CR(DEVD)₂ and Hoechst, and if a multi-wavelength filter option is

available on the fluorescence microscope, the dual staining properties can be examined. Hoechst Stain can be seen using a UV-filter with excitation at 365 nm and emission at 480 nm.

As AO exhibits a very broad emission range, one of several filter pairings may be used. The same excitation/emission pairing filters used to view CR(DEVD)₂ may be used: a 550 nm (540-560 nm) excitation and long pass >610 nm emission/barrier filter pairing. With this pairing, the lysosomes appear red. When illuminating with a blue light (480 nm) excitation filter, a green light (540-550 nm) emission/barrier filter combination works well. Lysosomes will appear yellowish green instead of red.

FIGURE 6: QUANTIFYING THE APOPTOTIC EFFECTS OF A DRUG AND CONTROLS ON PULMONARY ARTERY SMOOTH MUSCLE CELLS

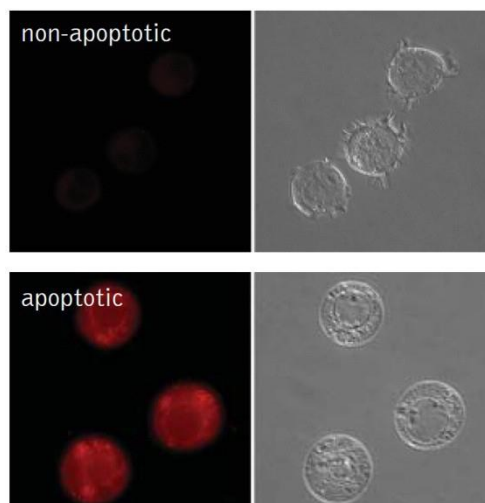


Enzo Life Sciences' CR(DEVD)₂ (BML-K1224), was used to quantify apoptosis via caspase 3&7 activity in human pulmonary artery smooth muscle cells (PASMC). Cells were treated with: a negative control condition (A); a drug that inhibits proliferation of PASMCs (B); or 1 mM H₂O₂ as a positive control to induce apoptosis (C). Cells were then labeled with CR(DEVD)₂ to detect caspases 3&7 and with Hoechst to stain nuclei blue. Cells in red (CR(DEVD)₂) contain active caspases 3&7 and have less intense blue nuclei (Hoechst) than healthy cells, which have bright blue nuclei. Only 11% of untreated cells (A) and 12% of drug-treated cells (B) were apoptotic (red) compared with 89% of cells treated with H₂O₂ (C). Although this drug inhibits proliferation of PASMCs, it does not induce apoptosis. Research done by Dr. Frederic Perros, et al., Universite Paris-Sud 11; INSERM U764, Clamart; INSERM U841, Hopital Henri-Mondor, Creteil, France¹⁶

8. Microscopy Analysis of Adherent Cells

1. Seed 104-105 cells onto a sterile coverslip in a 35 mm petri dish or onto chamber slides, or grow in a plate (Figure 2).
2. Grow cells until 80% confluent. This usually takes about 24 hours but will vary with each cell line.
3. Expose cells to the experimental conditions and create positive and negative controls (Section 2) or induce cells to undergo apoptosis (Section 3).
4. Reconstitute CR(DEVD)₂ to form the stock solution (Section 4):
 - a. Use 100 μ L DMSO to reconstitute the vial BML-KI224.
5. When ready to stain cells, dilute the stock 1:5 in diH₂O to form the CR(DEVD)₂ staining solution (Section 4):
 - a. Add 400 μ L diH₂O to dilute the vial BML-KI224.
 - b. Add CR(DEVD)₂ staining solution at approximately 1:30 and gently mix to ensure an even distribution of CR(DEVD)₂. For example, add 10 μ L staining solution to 300 μ L cells forming a final volume of 310 μ L. Do not add CR(DEVD)₂ to cells that will be stained with AO: add a placebo instead, such as diH₂O (Step 10).
7. Incubate 30-60 minutes at 37°C.
8. Remove the media from the cell monolayer surface and rinse twice with PBS, 1 minute per rinse. At this point, cells can be analyzed (Step 12) or stained with Hoechst (Step 9) or unstained cells can be labeled with AO (Step 10).

FIGURE 7: NEGATIVE VS. APOPTOTIC THP-1 CELLS



Using Enzo Life Sciences' CV-Caspase 3&7 Substrate Kit (BML-AK118-0001) to detect DEVDase activity in THP-1 cells, there is a clear differential between nonapoptotic (negative, top) and apoptotic (positive, bottom) cells. THP-1 cells were incubated with DMSO (top) or 1 μ M staurosporine (BML-EI156; bottom) for 3 hours at 37°C to induce DEVDase activity. Cell cultures were subsequently stained with CR(DEVD)₂ for 1 hour at 37°C. Left panels contain fluorescence images obtained using a Nikon Eclipse E800 photomicroscope equipped with a 100 W mercury lamp and excitation (510-560 nm) and emission (570-620 nm) filter pairings. Right panels contain the corresponding differential-interference-contrast (DIC) image (compare with Jurkat cells in Figure 8).

9. If cells are to be labeled with Hoechst Stain, add it at approximately 0.5% v/v. Add 1.55 μ L Hoechst to 310 μ L cells labeled with CR(DEVD)₂ and control samples. Incubate 5-10 minutes at 37°C. Go to Step 11.

10. Because of the overlap in emissions, dual staining of cells with both CR(DEVD)₂ and AO is not recommended; the dyes should be used separately. To stain cells that have not been exposed to CR(DEVD)₂:

a. Dilute AO at 1:2,000-1:200 (which is 0.05-0.5% v/v) into the final cell volume. For example, if using AO at 1.0 μ M in the final cell volume, it must be diluted 1:1,000. First dilute it 1:100 in PBS; e.g., add 10 μ L AO to 1,000 μ L PBS. Pipette the diluted AO to the cells at 1:10; e.g., add 35 μ L diluted AO to 315 μ L cell media.

b. Incubate 30 minutes at 37°C.

c. Remove the media from the cell monolayer surface. Rinse twice with PBS, 1 minute per rinse.

11. Mount the coverslip with cells facing down onto a drop of PBS. If a chamber-slide was used, pull off the plastic frame and add a drop of PBS to the cell surface and cover with a coverslip.

12. Observe cells using a fluorescence microscope equipped with an excitation filter of 550 nm (540- 560 nm) and a long pass >610 nm emission/barrier filter pair. Select a filter combination that best approximates these settings. Using these filters, positive cells stained with CR(DEVD)₂ will appear red with more brightly stained vacuoles and lysosomes.

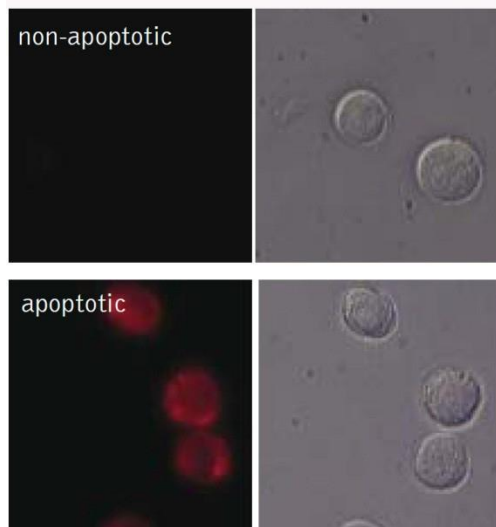
If samples were stained with both CR(DEVD)₂ and Hoechst, and if a multi-wavelength filter option is available on the fluorescence microscope, the dual

staining properties of the sample can be examined. Hoechst Stain can be seen using a UV-filter with excitation at 365 nm and emission at 480 nm.

As AO exhibits a very broad emission range, one of several filter pairs may be used. The same excitation/emission pairing filters used to view CR(DEVD)₂ may be used: an excitation filter of 550 nm (540-560 nm) and a long pass >610 nm emission/barrier filter. With this pairing, the lysosomes appear red.

When illuminating with a blue light (480 nm) excitation filter, a green light (540-550 nm) emission/barrier filter combination works well. Lysosomes will appear yellowish green instead of red.

FIGURE 8: DEVDASE ACTIVITY IN JURKAT CELLS



Following the same protocol outlined in Figure 7, Jurkat cells were analyzed to detect DEVDase activity. Negative cells do not fluoresce (top), while apoptotic cells fluoresce red using CR(DEVD)₂ (bottom). Left panels contain the fluorescence image; right panels contain the DIC image (compare with THP-1 cells in Figure 7).

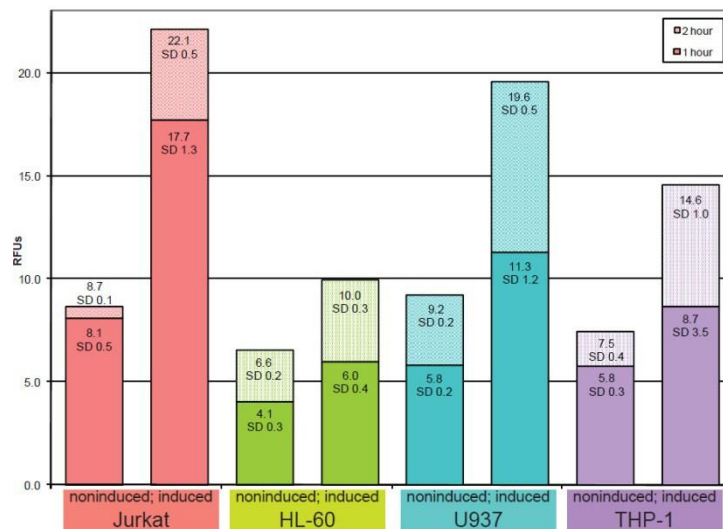
9. Fluorescence Plate Reader Staining Protocol

1. Prepare cell populations. Cell concentrations should be 2-8 x 10⁶ cells/mL. If this is too dense for the cell line, induce apoptosis first, then concentrate the cells and stain with CR(DEVD)₂. Cell concentration can be achieved by low speed centrifugation (<400 x g at RT) for 5 minutes.
2. Expose cells to the experimental conditions and create

positive and negative controls (Section 2) or induce cells to undergo apoptosis (Section 3).

3. Transfer 300 μL cell suspension into sterile tubes or a black microtiter plate. Do not use clear plates. Avoid bubbles. Larger cell volumes may also be used, but additional $\text{CR}(\text{DEVD})_2$ substrate will be required per sample.
4. When ready to label with the $\text{CR}(\text{DEVD})_2$ staining solution, cells should be at least 2×10^5 cells/100 μL aliquot (equal to 2×10^6 cells/mL) for each microtiter plate well.

FIGURE 9: QUANTIFICATION OF CASPASE ACTIVITY IN 4 CELL LINES USING A FLUORESCENCE PLATE READER



DEVDase activity in Jurkat, HL-60, U937, and THP-1 cells was quantified using Enzo Life Sciences' CV- Caspase 3&7 fluorogenic substrate, $\text{CR}(\text{DEVD})_2$ (BML- K1224), and analyzed with a fluorescence plate reader. Cells were incubated with 1 μM staurosporine (BML- E1156) or DMSO control for 3 hours (Jurkat cells) or 4 hours (HL-60, U937, THP-1 cells) at 37° C to induce caspase activity. Cells were incubated with $\text{CR}(\text{DEVD})_2$ for 1-2 hours to reveal the increase in intracellular DEVDase activity associated with apoptosis induction. DEVDase activity in each cell line is shown after 1 hour (solid bars) and 2 hour (spotted bars) exposures to the $\text{CR}(\text{DEVD})_2$. Cells were analyzed using a Molecular Devices Gemini XS fluorometric plate reader set at 590 nm excitation, 640 nm emission, with a 630 nm cut-off filter.

5. Reconstitute $\text{CR}(\text{DEVD})_2$ to form the stock solution (Section 4):
 - a. Use 100 μL DMSO to reconstitute the vial BML-K1224.
6. When ready to stain cells, dilute the stock 1:5 in diH_2O to form the $\text{CR}(\text{DEVD})_2$ staining solution (Section 4):

- a. Add 400 μL diH_2O to dilute the vial BML-KI224.
7. Add 20 μL $\text{CR}(\text{DEVD})_2$ staining solution directly to 300 μL cell sample. If different cell volumes are used, add $\text{CR}(\text{DEVD})_2$ staining solution at approximately 1:15. Due to sensitivity limitations, plate readers require a higher concentration of $\text{CR}(\text{DEVD})_2$ for detection compared with microscopes.
8. Gently mix the cells. This can be done by gently aspirating and expelling the cells with a pipette. To minimize cell shearing, cut the tip of the pipette to enlarge the hole.
9. Incubate cells for at least 60 minutes at 37°C protected from light. As cells settle to the bottom, gently resuspend them approximately every 20 minutes to ensure $\text{CR}(\text{DEVD})_2$ is evenly dispersed among all cells.
10. Read the 300 μL sample as one sample or split it into 3 wells of 100 μL each. If cells were stained in a tube, transfer 100-300 μL into a well of a black microtiter plate.
11. Measure the fluorescence intensity of the red fluorescent cresyl violet fluorophore. Set the plate reader to perform an endpoint read. $\text{CR}(\text{DEVD})_2$ has an optimal excitation and emission wavelength tandem of 592 nm and 628 nm, respectively. Select the filter pairings that best approximate these settings. If available, use a cut-off filter at 630 nm to filter out shorter wavelength excitation interference (Figure 9).

REFERENCES

1. Adamus, G., et al. Anti-enolase antibodies block enolase function and induce apoptosis of retinal cells. *Invest Ophthalmol Vis Sci.* 46:3478 (2005).
2. Boonacker, E., et al. Enzyme cytochemical techniques for metabolic mapping in living cells, with special reference to proteolysis. *J Histochem Cytochem.* 49:1473-1486 (2001).
3. Earnshaw, W. C., et al. Mammalian caspases: structural, activation, substrates, and functions during apoptosis. *Annu Rev Biochem.* 68:383-424 (1999).
4. Gordin, M., et al. C-Met and its ligand hepatocyte growth factor/ scatter factor regulate mature B cell survival in a pathway induced by CD74. *J Immunol.* 185:2020-2031 (2010).
5. Hwang, R. F., et al. Cancer-associated stromal fibroblasts promote pancreatic tumor progression. *Cancer Res.* 68:918- 926 (2008).
6. Idris, S. B., et al. Biocompatibility of polyester scaffolds with fibroblasts and osteoblast-like cells for bone tissue engineering. *J Bioactive Compatible Polymers.* 25:567-583 (2010).
7. Jang, S. W., et al. A selective TrkB agonist with potent neurotrophic activities by 7, 8-dihydroxyflavone. *PNAS.* 107: 2687-2692 (2010).
8. Johnson, D. E. Noncaspase proteases in apoptosis. *Leukemia* 14:1695-1703 (2000).
9. Khan, N., et al. Fisetin, a novel dietary flavonoid, causes apoptosis and cell cycle arrest in human prostate cancer LNCaP cells. *Carcinogenesis.* 29:1049-1056 (2008).
10. Kiang, J. G., et al. Inhibition of inducible nitric-oxide synthase protects human T cells from hypoxia-induced apoptosis. *Mol Pharmacol.* 73:738-747 (2008).
11. Lee, BW, et al. DEVDase detection in intact apoptotic cells using the cell permeant fluorogenic substrate, (z-DEVD)-2-cresyl violet. *BioTechniques.* 35(5):1080-1085 (2003).
12. Leist, M., et al. Triggering of apoptosis by cathepsins. *Cell Death Differ.* 8:324-326 (2001).
13. Lin, S., et al. Comparison of the toxicity of smoke from conventional and harm reduction cigarettes using human embryonic stem cells. *Toxicol Sci.* 118:202-212 (2010).
14. Morelli, M. B., et al. Characterization, expression, and functional activity of pituitary adenylate cyclase-activating polypeptide and its receptors in human granulosa-luteal cells. *J Clin Endocrinol Metab.* 93:4924-4932 (2008).
15. Nicholson, D. W. Caspase structure, proteolytic substrates, and function during apoptotic cell death. *Cell*

- Death Differ. 6:1028-1042 (1999).
16. Perros, F., et al. Platelet-derived growth factor expression and function in idiopathic pulmonary arterial hypertension. *Am J Respir Crit Care Med.* 178:81-88 (2008).
 17. Reed, J. C. Warner-Lambert/Parke-Davis award lecture: mechanisms of apoptosis. *Am J Pathol.* 157:1415-1430 (2000).
 18. Sanchez, C. G., et al. Activation of autophagy in mesenchymal stem cells provides tumor stromal support. *Carcinogenesis.* 32:964-972 (2011).
 19. Talanian, R. V., et al. Caspases as targets for anti-inflammatory and anti-apoptotic drug discovery. *J Med Chem.* 43:3351-3371 (2002).
 20. Thomas, A. J., et al. FOXD3 regulates the lineage switch between neural crest-derived glial cells and pigment cells by repressing MITF through a non-canonical mechanism. *Development.* 136:1849-1858 (2009).
 21. Thornberry, N. A. Caspases: key mediators of apoptosis. *Chem Biol.* 5:R97-R103 (1998).
 22. Thornberry, N. A., et al. A combinatorial approach defines specificities of members of the caspase family and granzyme. *B J Biol Chem.* 272:17907-17911 (1997).
 23. Van Noorden, C. J. F., et al. Ala-pro-cresyl violet, a synthetic fluorogenic substrate for the analysis of kinetic parameters of dipeptidyl peptidase IV (CD26) in individual living rat hepatocytes. *Anal Biochem.* 252:71-77 (1997).
 24. Voloboueva, L. A., et al. Mitochondrial protection attenuates inflammation-induced impairment of neurogenesis in vitro and in vivo. *J Neurosci.* 30:12242-12251 (2010).
 25. Wang, K. W. Calpain and caspase: can you tell the difference? *Trends Neurosci.* 23:20-26 (2000).
 26. Wei, Z., et al. Notch activation enhances the microglia-mediated inflammatory response associated with focal cerebral ischemia. *Stroke.* 42:2589-2594 (2011).



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:

www.enzolifesciences.com