

FAM-FLICA® Caspase Assay Kits



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

1. Introduction

FLICA® is a powerful method to assess cell death by detecting apoptosis *in vitro*. ImmunoChemistry Technologies' (ICT) FLICA® probes are non-cytotoxic Fluorescent Labeled Inhibitors of CAspases that covalently bind to active caspase enzymes^{1,2}. FLICA® measures the intracellular process of apoptosis instead of a side-effect, such as the turn-over of phosphatidyl serine, and eliminates the incidence of false positives that often plagues both Annexin V and TUNEL staining.

To use FLICA®, add it directly to the cell culture media, incubate, and wash. FLICA® is cell-permeant and will efficiently diffuse in and out of all cells. If there is an active caspase enzyme inside the cell, it will covalently bind with FLICA® and retain the green fluorescent signal within the cell. Unbound FLICA® will diffuse out of the cell during the wash step. Apoptotic cells will retain a higher concentration of FLICA® and fluoresce brighter than non-apoptotic cells. There is no interference from pro-caspases or inactive forms of the enzymes. If the treatment is causing cell death via apoptosis, apoptotic cells will have an elevated level of caspase activity relative to non-apoptotic cells or negative control cells and fluoresce with FLICA®.

Apoptosis is an evolutionarily conserved process of programmed cell suicide. It is centered on a cascade of proteolytic enzymes called caspases that are triggered in response to pro-apoptotic signals. Once activated, caspases cleave protein substrates leading to the eventual disassembly of the cell²⁵. Caspases have been identified in organisms ranging from *C. elegans* to humans. Mammalian caspases play distinct roles in both apoptosis and inflammation. In apoptosis, effector caspases (-3, -6, and -7) are responsible for proteolytic cleavages that lead to cell disassembly. Initiator caspases (-8, -9, and -10) regulate apoptosis upstream. Caspase-1 is associated with inflammasome activity and takes on the role of a key housekeeping enzyme in its conversion of pro-IL-1 β protein into the active IL-1 β cytokine. (Use FLICA® kits #98 and 9122 to detect caspase-1.)

Like the majority of other proteases, caspases are synthesized as pro-form precursors that undergo proteolytic maturation, either autocatalytically or in a cascade by enzymes with similar specificity¹⁵. Active caspase enzymes consist of two large (~20 kD) and two small (~10 kD) subunits that non-covalently associate to form a two heterodimer, tetrameric active caspase^{24,29,31}.

Activated caspase enzymes cleave proteins by recognizing a 3 or 4 amino acid sequence that must include an aspartic acid (D)

residue in the P1 position. This C-terminal residue is the target for the cleavage reaction at the carbonyl end²⁸. Each FLICA® probe contains a 3 or 4 amino acid sequence that is targeted by different activated caspases. This target sequence is sandwiched between a green fluorescent label, carboxyfluorescein (FAM), and a fluoromethyl ketone (FMK). A caspase enzyme cannot cleave the FLICA® inhibitor probe; instead, it forms an irreversible covalent bond with the FMK on the target sequence and becomes inhibited from further enzymatic activity. ICT's poly caspase FLICA® probe, FAM-VAD-FMK, may be used as a general reagent to detect apoptosis as it is recognized by all types of activated caspases. To more specifically target a particular caspase enzyme, use one of ICT's specialized FLICA® reagents. These reagents contain different amino acid target sequences preferred by each caspase: caspase-1 (YVAD), -2 (VDVAD), -3/7 (DEVD), -6 (VEID), -8 (LETD), -9 (LEHD), -10 (AEVD), and -13 (LEED). Some of these reagents are also available with a red or far red fluorescent label.

FLICA® has been known to detect caspase activity in human, rabbit, rat, mice, drosophila, squid, paramecium, and yeast cell lines, among others. FLICA® can be used to label suspension or adherent cells, thin tissue sections, and frozen sections. After labeling with FAM-FLICA®, cells can be fixed or frozen. For tissues that will be paraffin-embedded after labeling, use ICT's red sulforhodamine SR-FLICA® probes; do not use the green FAM-FLICA® probes as the green FAM dye will be quenched during the paraffin embedding process.

Cells labeled with FAM-FLICA® can be counter-stained with reagents such as the red vital stains Propidium Iodide (included in FAM-FLICA® kits) and 7-AAD (catalog # 6163) to distinguish apoptosis from necrosis. Nuclear morphology may be concurrently observed using Hoechst 33342, a blue DNA binding dye (included in FLICA® kits). Cells can be viewed directly through a fluorescence microscope (Figures 1-3, 9, and 10), or the fluorescence intensity can be quantified using a flow cytometer (Figures 4-7) or fluorescence plate reader (Figure 8). FAM-FLICA® optimally excites at 488-492 nm and has a peak emission at 515-535 nm. FLICA® is for research use only. Not for use in diagnostic procedures.

*Learn more about all of ICT's products at
www.immunochemistry.com or call 1-800-829-3194.*

2. FAM-FLICA Kits:

Caspase	Inhibitor Reagent (part#)	Trial	Standard
Poly Caspase	FAM-VAD-FMK (637)	91	92
Caspase-1	FAM-YVAD-FMK (655)	97	98
Caspase-2	FAM-VDVAD-FMK (680)	918	919
Caspase-3/7	FAM-DEVD-FMK (653)	93	94
Caspase-6	FAM-VEID-FMK (654)	95	96
Caspase-8	FAM-LETD-FMK (656)	99	910
Caspase-9	FAM-LEHD-FMK (677)	912	913
Caspase-10	FAM-AEVD-FMK (682)	922	923
Caspase-13	FAM-LEED-FMK (683)	929	930

3. Kit Contents

Trial size kits contain:

- 1 vial of FAM-FLICA caspase inhibitor reagent
- 1 bottle of 10X Apoptosis Wash Buffer (15 mL) #635
- 1 bottle of Fixative (6 mL) #636
- 1 vial of Propidium Iodide, 250 µg/mL (1 mL) #638
- 1 vial of Hoechst 33342, 200 µg/mL (1 mL) #639

Standard size kits contain:

- 4 vials of FAM-FLICA caspase inhibitor reagent
- 1 bottle of 10X Apoptosis Wash Buffer (60 mL) #634
- 1 bottle of Fixative (6 mL) #636
- 1 vial of Propidium Iodide, 250 µg/mL (1 mL) #638
- 1 vial of Hoechst 33342, 200 µg/mL (1 mL) #639

4. Storage

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

5. Safety Data Sheets (SDS)

Safety data sheets are available online at www.immunochemistry.com or by calling 1-800-829-3194 or 952-888-8788.

6. Recommended Materials

- DMSO, 50 µL per vial to reconstitute FLICA
- DiH₂O, 135-540 mL to dilute 10X Apoptosis Wash Buffer
- Phosphate buffered saline (PBS) pH 7.4, up to 100 mL, to dilute FLICA and handle cells
- FBS and/or BSA to add to the buffer when handling cells
- Cultured cells or tissues treated with the experimental conditions ready to be labeled
- Reagents to induce caspase activity and create controls, such as staurosporine (catalog #6212) or camptothecin (catalog #6210)
- 90% ETOH or 3% formaldehyde to create live/dead controls for Propidium Iodide staining
- Black 96-well microtiter plate, flat bottom, non-treated, non-sterile (ICT catalog #266). If using a bottom reading instrument, use a plate with black walls and a clear bottom. If culturing cells in the plate, use a sterile black tissue culture plate.
- Hemocytometer
- Centrifuge at <400 g
- 15 mL polypropylene centrifuge tubes (1 per sample)

7. Detection Equipment

The assay can be analyzed with a:

- Fluorescence microscope
- Fluorescence plate reader
- Flow cytometer

Use filter pairings that best approximate these settings:

- FAM-FLICA optimally excites at 488-492 nm and has a peak emission at 515-535 nm.
- View Propidium Iodide (PI) under a long pass filter with the excitation at 488-492 nm, emission >610 nm; nuclei-bound PI has a maximum emission at 617 nm (Section 13).
- Hoechst 33342 can be seen using a UV-filter with excitation at 365 nm and emission at 480 nm (Section 14).

8. Experimental Preparation

Staining apoptotic cells with FLICA can be completed within a few hours. However, FLICA 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 treatment or apoptosis induction process, which typically requires a 2-6 hour incubation at 37°C based on the cell line and concentration. Create cell populations, such as:

- a. Cells that were exposed to the experimental treatment
- b. A negative control population of cells that received a placebo treatment

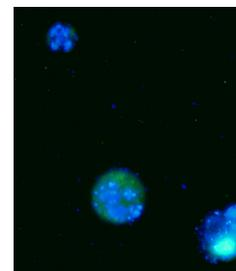
As FLICA detects the presence of catalytically active forms of caspase enzymes, plan the experiment so that FLICA will be diluted and administered at the time when caspases are expected to be activated in the cells.

The recommended volume of 30X FLICA is 10 µL per 300 µL of cells at 5 x 10⁵ cells/mL, but the amount may vary based on the experimental conditions and the instrument used for analysis. Each investigator should adjust the amount of FLICA to accommodate the particular cell line and research conditions.

Culture cells to a density optimal for the specific experimental conditions or apoptosis induction protocol. Cell density should not exceed 10⁶ cells/mL. Cells cultivated in excess of this concentration may begin to naturally enter apoptosis. An initial experiment may be necessary to determine when and how much FLICA to use as the resulting positive signal is a direct measurement of caspase activity occurring during the incubation period.

Figure 1: Microscopy Analysis of THP-1 Suspension Cells

Human monocytic leukemia THP-1 cells were dually stained with ICT's green FAM-FLICA poly caspase inhibitor reagent, FAM-VAD-FMK (kit #92), and a blue DNA stain, Hoechst 33342. Cells were incubated with 1 µM staurosporine for 3 hours at 37°C to induce apoptosis. Cells were then labeled with FAM-VAD-FMK for 60 minutes at 37°C. Cells were washed, then Hoechst stain was added and incubated for 5 minutes. Wet-mount slides were prepared and two photos were taken and superimposed. Caspase activity (green) was detected using a band pass filter (excitation at 488 nm; emission at 520 nm). Nuclear staining by Hoechst 33342 was revealed using a UV-filter (excitation at 365 nm, emission at 480 nm). Only one cell of the three cells is apoptotic (middle) – it is stained positively for caspase activity with FLICA FAM-VAD-FMK. It also has many bright blue spots from the Hoechst stain, indicating that the DNA is breaking down and the cell is beginning to die. The lack of green staining and the concentrated blue DNA in the lower right cell indicate it is alive (not apoptotic). The upper left cell is necrotic (no green, scattered blue). Data courtesy of Dr. Brian W. Lee, ICT.



9. Controls

Create experimental samples and control cell populations:

- Treated experimental population(s): cells exposed to the experimental condition(s).
- Negative control: non-treated cells grown in a normal cell culture environment.
- Positive control: cells induced to undergo apoptosis using a known caspase activation protocol.

The induced positive cell population and negative control cell population tubes should come from a common pool of cells and contain similar quantities of cells. Create negative controls by culturing an equal volume of non-induced cells for every labeling condition. For example, if labeling with FLICA and Hoechst stain, make 8 populations:

- 1&2. Unlabeled: induced and non-induced
- 3&4. FLICA-labeled: induced and non-induced
- 5&6. FLICA-labeled and Hoechst-labeled: induced and non-induced
- 7&8. Hoechst-labeled: induced and non-induced

If analyzing cells with a flow cytometer, create Propidium Iodide (PI) instrument controls using formaldehyde or ETOH to compensate for bleed-over of the PI signal into FL-1 (Sections 13 and 19).

10. Apoptosis Induction

Prior to commencing the experiment, determine a reproducible method for obtaining a positive control by triggering caspase activity. This process varies significantly with each cell line. For example, apoptosis may be induced with 2-4 $\mu\text{g}/\text{mL}$ camptothecin or 1-2 μM staurosporine for >4 hours.

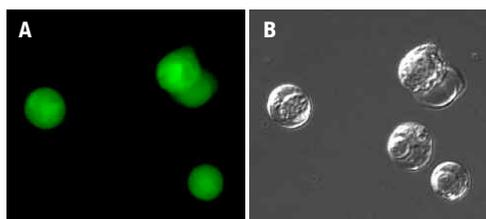
11. Preparation of FLICA

FLICA is supplied as a lyophilized powder that may be slightly visible as an iridescent sheen inside the vial. Protect from light and use gloves when handling. Because the 30X FLICA solution must be used immediately, prepare it just before staining.

- Reconstitute each vial of FLICA with 50 μL DMSO to form the 150X stock. The stock solution should be colorless to light yellow. Once reconstituted, it may be stored at $\leq -20^\circ\text{C}$ for 6 months protected from light and thawed no more than twice during that time.
- Immediately prior to addition to the samples and controls, dilute FLICA 1:5 by adding 200 μL PBS to each vial to form the 30X FLICA solution. Use 30X FLICA within 30 minutes of dilution into aqueous buffers.

Figure 2: Caspase Activity in Jurkat Cells

Jurkat cells (T lymphocytes) were labeled with ICT's green FAM-FLICA poly caspase inhibitor FAM-VAD-FMK (kit #92) and viewed under a fluorescence microscope. The grey DIC image (B) reveals five cells in the field of view, but only four of them fluoresce green (A). Four out of five cells are apoptotic and have active caspases present (green). Data courtesy of Dr. Brian W. Lee, ICT.



12. Preparation of 1X Apoptosis Wash Buffer

ICT's Apoptosis Wash Buffer (catalog #634 and #635, AWB) is an isotonic solution used to wash cells following exposure to FLICA. It contains mammalian proteins to stabilize cells stained with FLICA and sodium azide to retard bacterial growth (1X Apoptosis Wash Buffer contains 0.01% w/v sodium azide). Cell culture media containing FBS and other additives may be used to wash cells instead of Apoptosis Wash Buffer.

- 10X Apoptosis Wash Buffer may form precipitates during cold storage. If this happens, gently warm it until all crystals have dissolved. Do not boil.
 - Dilute 10X Apoptosis Wash Buffer 1:10 in diH₂O. For example, add 15 mL 10X Apoptosis Wash Buffer to 135 mL diH₂O for a total of 150 mL.
- 1X Apoptosis Wash Buffer may be stored at 2-8°C and used within 1 week or frozen and used within 6 months.

13. Propidium Iodide

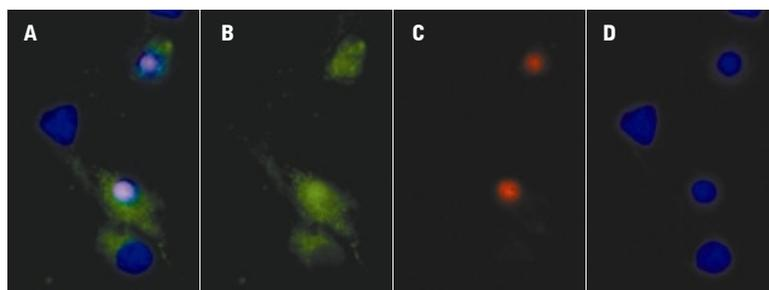
Propidium Iodide (PI, catalog #638) is used to distinguish between living and dead cells by staining necrotic, dead, and membrane-compromised cells red. PI is an intercalating fluorescent reagent that stoichiometrically binds every four to five base pairs of DNA. PI is membrane impermeant, which prevents it from reaching the DNA in viable cells, thus allowing the identification of dead cells in a population.

Upon binding to DNA, the fluorescence intensity potential of PI is enhanced 20-30 fold. Nucleic acid-bound PI has a red-shifted absorbance/excitation maximum of 535 nm and an emission maximum of 617 nm. PI efficiently excites at 488-492 nm. Its excitation and emission spectra allow for efficient analysis using fluorescence microscopy or flow cytometry. When reading with FAM in FL-1, PI may be read in FL-2 or FL-3 (it may be easier to compensate using FL-3 as there is less bleed over). PI is provided in a soluble, ready-to-use formulation (1 mL at 250 $\mu\text{g}/\text{mL}$).

- Warning:** Propidium Iodide is a mutagen. It may cause serious eye irritation. Gloves, protective clothing, and eye wear are strongly recommended. When disposing, flush sink with copious amounts of water; see SDS for further information.

Figure 3: Cell Death in Primary Rat Hippocampal Neurons

ICT's FAM-FLICA caspase-3/7 inhibitor reagent, FAM-DEVD-FMK (kit #94), was used to assess cell death in primary rat hippocampal neurons. Subjects were first-generation descendants of Sprague-Dawley albino rats. Hippocampi from PND 0 male pups were used for primary cultures of hippocampal neurons. Cells were plated on 25-mm poly-L-lysine-coated coverslips at 300,000 cells per coverslip, and cells were used at 4 or 8 days *in vitro*. Image A is a composite of FLICA (B), PI (C), and Hoechst (D). In A, 4 cells are revealed by labeling DNA with Hoechst (blue, D), 3 out of the 4 cells are caspase-positive (green, B), and 2 of those cells are also membrane-compromised (red, C). 3 of the cells fluoresce green with FLICA FAM-DEVD-FMK (B); they are caspase-positive and undergoing apoptosis. 1 of the cells is in early apoptosis as it is green but not red. 2 of the FLICA-positive cells are also PI-positive (red, C) meaning they are becoming membrane compromised and are in the late stages of apoptosis rather than necrosis. Data courtesy of Dr. Z. Kahraman Akozer, University of Maryland.



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

PI can be used with FAM-FLICA to identify four populations of cells: living; early apoptotic; late apoptotic; and necrotic (Figures 3, 5, 6). If analyzing PI-labeled cells using flow cytometry (Section 19), create PI instrument controls using formaldehyde or ETOH to kill cells. An apoptosis-induction method (Section 10) is not as effective as a solvent treatment at creating PI-positive cell controls for instrument calibration because apoptosis-induced samples may not have enough late-stage apoptotic cells that have become membrane-compromised to stain positive for PI. To set up PI instrument controls:

- Label 2 centrifuge tubes:
 - PI-negative (live)
 - PI-positive (dead)
- Add 1-5 x 10⁵ non-induced, healthy cells to each tube.
- Centrifuge at 200 x g for 5 minutes at room temperature (RT) to pellet cells; remove supernatants.
- To create the PI-negative control (live cells), resuspend the cells in 300 μ L 1X AWB or PBS+1% BSA to maintain the integrity of the cell membrane.
- To create the PI-positive control (membrane-compromised dying cells):
 - Use formaldehyde: resuspend cells in 300 μ L 3% v/v formaldehyde (in 97% PBS or 1X AWB). Incubate 30 minutes on ice and then wash cells (Step 6).
 - Or use ETOH: resuspend cells in 300 μ L 90% ETOH (in 10% PBS or 1X AWB). Gently vortex for 30 seconds then wash cells (Step 6).
- Add 1 mL 1X AWB or PBS+1% BSA.
- Centrifuge at 200 x g for 5 minutes; remove supernatants.
- Resuspend in 500 μ L 1X AWB or PBS+1% BSA.
- If ETOH was used, add 500 μ L non-induced, healthy cells to the tube of ETOH-killed cells and mix; use roughly the same number of cells as originally included in Step 2. This will create a sample with distinct positive and negative peaks.
- Add PI at 0.5% v/v. For example, add 2.5 μ L PI to a 500 μ L sample. Incubate 5 minutes, protected from light at RT.
- Read immediately on the flow cytometer and compensate bleed-over of the red PI signal from FL-2 or FL-3 into FL-1. It may be easier to compensate PI spillover into FL-1 when read in FL-3.

14. Hoechst 33342

Hoechst 33342 is a cell-permeant nuclear stain that emits blue fluorescence when bound to double stranded DNA. It is used to stain the nuclei of living or fixed cells, to distinguish condensed pyknotic nuclei in apoptotic cells, and for cell cycle studies. Hoechst 33342 is provided ready-to-use at 200 μ g/mL. Hoechst 33342 can be used with FAM-FLICA and PI to label the nuclei of live, dying, and apoptotic cells.

When bound to nucleic acids, it has a maximum absorbance at 350 nm and a maximum emission at 480 nm. It is revealed under a microscope using a UV-filter with excitation at 365 nm and emission at 480 nm.

- Warning:** Hoechst 33342 is a potential mutagen. It may be irritating to respiratory system and skin. Gloves, protective clothing, and eye wear are strongly recommended. When disposing, flush sink with copious amounts of water. See SDS for further information.

15. Fixative

ICT's Fixative (catalog #636) is a formaldehyde solution designed to cross-link and aggregate intracellular components. If the stained cell populations cannot be evaluated immediately after labeling with FLICA, add Fixative at a ratio of 1:10. For example, add 100 μ L Fixative to 900 μ L cells. Never add Fixative until all the staining and final wash steps have been completed. Fixed cells may be stored on ice or at 4°C for up to 24 hours, protected from light.

ICT's Fixative will not interfere with the carboxyfluorescein (FAM) label. Do not use absolute ethanol- or methanol-based fixatives as they will inactivate the FAM-FLICA label. Do not fix cells that will be stained later with Propidium Iodide or Hoechst 33342.

- Warning:** Fixative is toxic: danger exists of very serious irreversible effects through inhalation, by contact with skin, or if swallowed. Gloves, protective clothing, and eye wear are strongly recommended. When disposing, flush sink with copious amounts of water; see SDS for further information.

16. Staining Protocol for Suspension Cells

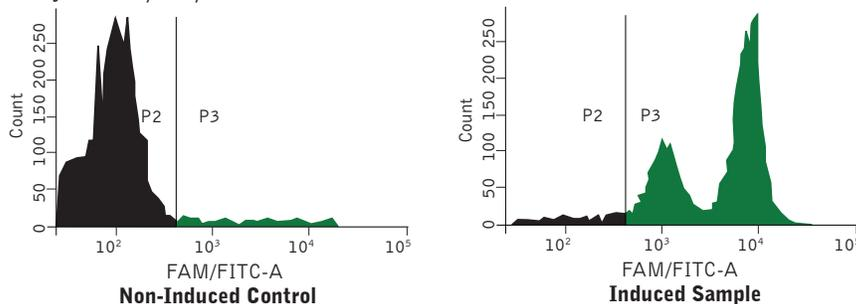
Prepare experimental and control cell populations. Ideally, cell concentration should be 3-5 x 10⁵ cells/mL. The concentration should not exceed 10⁶ cells/mL, as cells cultivated in excess of this concentration may begin to naturally enter apoptosis. Just prior to staining with FLICA, cells may need to be concentrated to 2-5 x 10⁶ cells/mL as both microscopy and plate reader analysis methods (Sections 17 and 20) require high cell concentrations. Start with a larger volume of cells at 3-5 x 10⁵ cells/mL (which is a typical density for cell culture) and then concentrate cells and resuspend to 300 μ L per sample when ready for FLICA staining.

- Expose cells to the experimental or control condition. If analyzing with a flow cytometer, set aside four populations to create instrument controls with PI-positive and PI-negative cells, and FLICA-induced and FLICA-non-induced cells
- If analyzing with a fluorescence microscope or plate reader, concentrate cells to 2-5 x 10⁶ cells/mL just prior to FLICA staining. Fluorescence microscopy requires an excess of 2 x 10⁶ cells/mL to obtain 5 - 20 cells per image field. Flow cytometry can analyze samples at 3-5 x 10⁵ cells/mL.
- Transfer 290 μ L cells into fresh tubes.

Figure 4: Single Color Analysis via Flow Cytometry

Jurkat cells were treated with DMSO, a negative control (left), or staurosporine, an apoptosis-inducing agent (right), for 4 hours, then stained with ICT's green poly caspase inhibitor probe, FAM-VAD-FMK (kit #92), for 1 hour. Cells were washed twice and read on a flow cytometer. The negative control exhibited caspase activity in only 3.3% of the cell population (P3, left histogram), whereas treatment with staurosporine induced caspase activity in 94.8% of the experimental cells (P3, right histogram). This is a ratio of 28:1. Data courtesy of Ms. Tracy Hanson, ICT, 10G5.

% of Cells	P2 Negative	P3 Positive
Negative, Non-Induced (left)	87.5%	3.3%
Positive, Induced (right)	4.8%	94.8%
Ratio		28:1



4. Add 10 μL 30X FLICA solution, forming a final volume of 300 μL . If different cell volumes were used, add 30X FLICA at a ratio of 1:30. Mix by gently flicking the tubes. The amount of FLICA should be optimized for each cell line and experimental condition.
5. Incubate cells at 37°C protected from light. The incubation period may range from 30 minutes to several hours and should be optimized for each cell line and experimental condition. As cells may settle on the bottom of the tubes, gently resuspend them by swirling cells every 20 minutes to ensure an even distribution of FLICA.
6. If cells are to be analyzed with a microscope, cells may be dually stained with Hoechst. Add Hoechst at 0.5% v/v and incubate 5 minutes at 37°C. For example, if the cell suspension is at 300 μL , add 1.5 μL Hoechst.
7. Add 2 mL 1X Apoptosis Wash Buffer and gently mix.
8. Centrifuge at 200 x g for 5 minutes at RT.
9. Carefully remove and discard supernatants. Gently vortex the pellets to disrupt clumping. Resuspend in 1 mL 1X Apoptosis Wash Buffer and gently mix.
10. Centrifuge cells at 200 x g for 5 minutes at RT.
11. Carefully remove and discard supernatants. Gently vortex pellets to disrupt clumping.
 - If analyzing with a fluorescence microscope, go to Section 17.
 - If using a flow cytometer, go to Section 18 or 19.
 - If using a plate reader, go to Section 20.

17. Microscopy Analysis of Suspension Cells

Follow Section 16, Steps 1-11. See Figures 1-2.

12. Resuspend cells in 300-500 μL 1X Apoptosis Wash Buffer and place on ice. At this point, the cells may be stained with Propidium Iodide (PI) for bicolor analysis (Step 13), fixed for future viewing (Step 14), or observed immediately (Step 15).
13. To identify dead cells by staining with PI, add 1.5 μL PI to the 300 μL cell suspension. If different volumes were used, add it at 0.5% v/v. Incubate 5 minutes at 37°C.
 - a. Wash cells to remove excess PI from the media. Centrifuge at 200 x g for 5 minutes at RT.
 - b. Carefully remove and discard supernatants. Gently vortex pellets to disrupt clumping.
 - c. Resuspend cells in 300 μL 1X Apoptosis Wash Buffer and gently mix. Go to Step 14 or 15.
14. If not viewing immediately, cells may be fixed for viewing up to 24 hours later. Add 30 μL Fixative. If cells were resuspended in a different volume, add it at a ratio of 1:10.
 - a. Incubate 15 minutes at RT in the dark.
 - b. Place cells onto a microscope slide and allow to dry.
 - c. Briefly wash cells with PBS.
 - d. Cover cells with mounting media and coverslip.
 - e. Store slides at 2-8°C for up to 24 hours.
15. To view cells immediately, place 1 drop of cell suspension onto a microscope slide and cover with a coverslip.
16. Observe cells under a fluorescence microscope using a bandpass filter (excitation 490 nm, emission >520 nm) to view green fluorescence. Cells bearing active caspase enzymes covalently coupled to FLICA appear green (Figures 1-3, 6, and 9). Hoechst stain can be seen using a UV-filter with excitation at 365 nm and emission at 480 nm (Figures 1 and 3). View nucleic-acid-bound PI under a long pass filter with the excitation at 490 nm, emission >610 nm.; Nuclei-bound PI has a maximum emission at 617 nm (Figures 3 and 6).

Figure 5: Bicolor Staining of Rabbit Cells

ICT's poly caspase inhibitor FLICA reagent, FAM-VAD-FMK (kit #92), was used to assess apoptosis in rabbit cells (both suspension and adherent cells). Cells were grown to 3×10^6 cells per sample and treated with a placebo (top) or a condition that induced apoptosis (bottom). Cells were stained with FAM-VAD-FMK, washed, stained with the red vital stain Propidium Iodide (PI), and analyzed using two-color flow cytometry. Dot plots were set up to detect caspase activity (green, FL-1) on the X-axis and necrosis (red, FL-2) on the Y-axis. Four populations of cells were detected: (A) unstained live cells do not fluoresce; (D) cells in early apoptosis fluoresce green with FAM-FLICA; (C) cells in late apoptosis are dually stained with FAM-FLICA and PI: they fluoresce green (active caspases) and red (the cell membrane has been permeabilized); and (B) necrotic cells fluoresce red. Cells became necrotic when treated with the placebo (top, B and D), while many cells entered early and late apoptosis when treated with the inducer (bottom, C and D). Compare data with Figure 6. Data courtesy of JH at HMC.

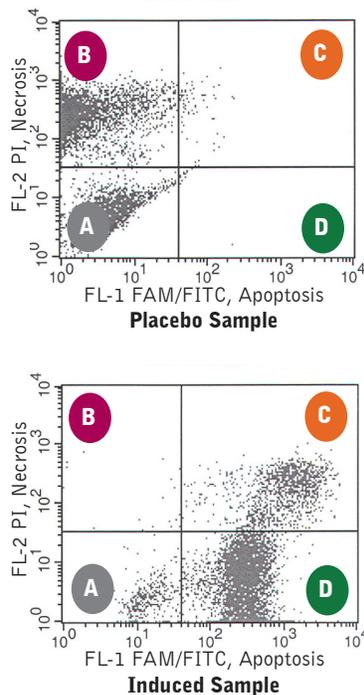
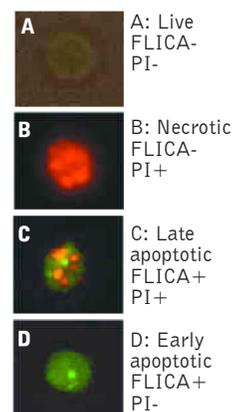
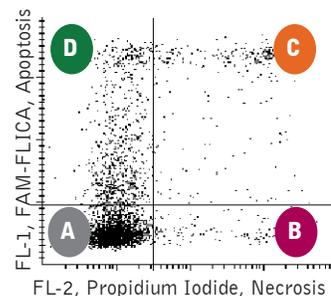


Figure 6: Four Populations of HL-60 Cells

HL-60 cells (human promyelocytic leukemia) were treated with a drug, then stained with ICT's poly caspase inhibitor reagent, FLICA FAM-VAD-FMK (kit #92), and a red vital stain, Propidium Iodide (PI). Cells were analyzed on a scanning laser cytometer to detect red on the X-axis (necrosis) and green on the Y-axis (apoptosis). Four populations of cells were detected: (A) unstained live cells do not fluoresce; (D) cells in early apoptosis fluoresce green with FAM-FLICA; (C) cells in late apoptosis are dually stained with FAM-FLICA and PI: they fluoresce green (active caspases) and red (the cell membrane is compromised); and (B) necrotic cells fluoresce red. In this experiment, the drug triggered the caspase cascade rather than being overtly toxic to the cells. Compare data with Figure 5. Data courtesy of Dr. Zbigniew Darzynkiewicz, Brander Cancer Center, NY.



FOR RESEARCH USE ONLY.

NOT FOR USE IN DIAGNOSTIC PROCEDURES.

18. Single Color Flow Cytometry Analysis of Suspension Cells

Follow Section 16, Steps 1-5 and 7-11 (omit Hoechst staining in Step 16.6). See Figure 4.

- Resuspend cells in 300 μL 1X Apoptosis Wash Buffer and place on ice.
- For single-color analysis, use a 15 mW argon ion laser at 488 nm. Measure fluorescein on the FL-1 channel. Generate a histogram with the log FL-1 on the X-axis versus the number of cells on the Y-axis. Caspase negative (FLICA-) cells will occur in the lower log fluorescence output decades of the FL-1 (X) axis, whereas caspase-positive (FLICA+) cells will appear as a shoulder on the right side (brighter) or separate peak on the right side of the negative peak histogram (Figure 4).

19. Multicolor Flow Cytometry Analysis of Suspension Cells

Follow Section 16, Steps 1-5 and 7-11 (omit Hoechst staining, Step 16.6). See Figures 5-7.

To address compensation issues and set up the flow cytometer, prepare two sets of instrument controls: live and killed cells that are only stained with PI (Section 13); and induced and non-induced cells that are only stained with FLICA. These controls are needed to adjust the instrument PMT's to separate PI-positive and PI-negative samples and to compensate for bleed-over of the red PI signal from FL-2 or FL-3 into FL-1. They will also help to clearly differentiate the FLICA-positive population from the FLICA-negative population and compensate bleed-over of the green FAM-FLICA signal from FL-1 into FL-2 or FL-3. If using an instrument that does not have the option to adjust PMT's, like the Accuri C6, these control populations can help set the positive gate for the PI-positive and FLICA-positive cells prior to reading dual-labeled samples.

While setting up the PI controls (Section 13), continue working with the test samples and experimental controls as prepared in Section 16.

- Resuspend cells in 400 μL 1X Apoptosis Wash Buffer.
- Set aside a control of FLICA-stained induced cells that does not contain PI. FLICA-stained cells may be stored at 2-8°C for up to 4 hours protected from light, depending upon the cell line.
- Stain cells with 2 μL PI, mix, and put on ice. Analyze as quickly as possible (within 30 minutes) as PI-stained cells may begin to have toxicity issues. If not analyzing PI-stained cells within 30 minutes, wash and fix:
 - Wash cells to remove the PI from the media to prevent false-positives as any excess PI will go into all cells after fixation. Centrifuge at 200 x g for 5 minutes at RT.
 - Carefully remove and discard supernatants. Gently vortex the pellets to disrupt clumping.
 - Resuspend cells in 300 μL 1X Apoptosis Wash Buffer and gently mix.
 - Add 30 μL Fixative. If cells were resuspended in a different volume, add Fixative at a ratio of 1:10.
 - Incubate 15 minutes at RT in the dark. Store fixed cells at 2-8°C for up to 24 hours protected from light.

- Set up the instrument compensation.
 - Read the PI-positive and PI-negative controls (Section 13) to compensate bleed-over of the red PI signal from FL-2 or FL-3 into FL-1.
 - Read the FLICA-only-positive and FLICA-only-negative controls to compensate bleed-over of the green FAM-FLICA signal from FL-1 into FL-2 or FL-3.
- When ready to read the samples for bicolor analysis, measure carboxyfluorescein (FAM) on the FL-1 channel and red fluorescence (PI) on the FL-2 or FL-3 channel. Generate a log FL-1 versus log FL-2 or FL-3 dot plot (Figures 5 and 6). This will reveal four populations of cells:
 - Cells in early apoptosis fluoresce green with FAM-FLICA.
 - Cells in late apoptosis are dually stained with FAM-FLICA and PI; they fluoresce green (they have active caspases) and red (the cell membrane has permeabilized).
 - Necrotic cells fluoresce red.
 - Unstained live cells do not fluoresce.

20. Fluorescence Plate Reader Analysis of Suspension Cells

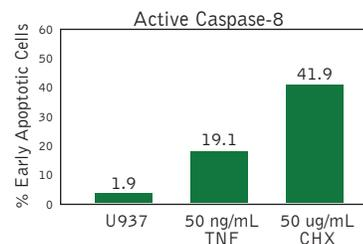
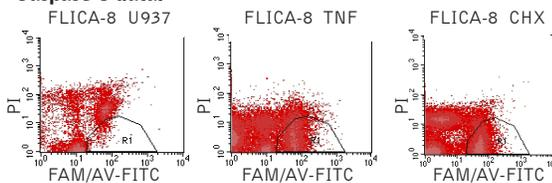
Follow Section 16, Steps 1-5 and 7-11 (omit Hoechst staining, Step 16.6). See Figure 8.

- Resuspend cells in 500 μL PBS.
- Determine the concentration and compare the cell density of each sample. The non-induced population may have more cells than the induced population, as some apoptotic cells in the induced samples may be lost during the wash steps. Adjust the volume of the cell suspensions to equalize the cell density. When ready to read, cells should be $>3 \times 10^6$ cells/mL.
- Pipette 100 μL per well into a black microtiter plate. Analyze at least 2 aliquots per sample. Do not use clear plates. Avoid bubbles.
- Perform an endpoint read. Set the excitation wavelength at 488 nm and the emission wavelength to 530 nm; use a cut-off filter of 515 nm. FAM-FLICA has an optimal excitation range of 488 - 492 nm and emission range of 515 - 535 nm (Figure 8).

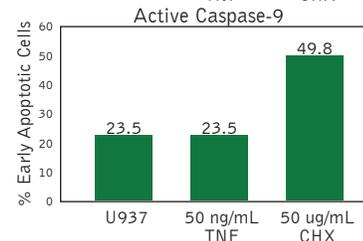
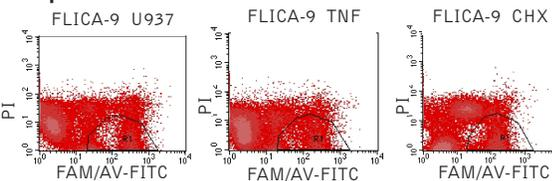
Figure 7: Flow Cytometry Analysis

ICT's caspase-8 and caspase-9 FAM-FLICA kits were used to detect early apoptosis in U937 cells, a histiocytic lymphoma cell line, upon treatment with TNF or cycloheximide for 24 hours. 200 μL of U937 cells at 10^6 cells/mL were pipetted into a 96-well U-bottom plate. Apoptosis was induced by treating cells with either 50 ng/mL TNF or 50 $\mu\text{g}/\text{mL}$ cycloheximide (CHX) for 24 hours. Cells were labeled by adding 6 μL 30X FLICA caspase-8 inhibitor reagent (FAM-LETD-FMK, kit #910) or FLICA caspase-9 inhibitor reagent (FAM-LEHD-FMK, kit #913) for 1 hour. Cells were washed three times, stained with propidium iodide (PI), and analyzed with a BD FACS. Caspase-9 was more active in these cells than caspase-8. Very few of the cells became necrotic (PI-positive). Data courtesy of JM at SRI.

Caspase-8 data:



Caspase-9 data:



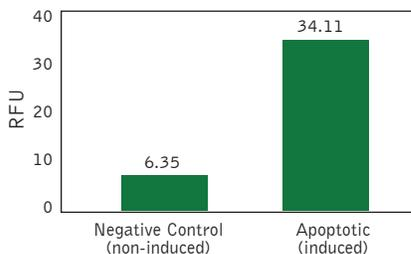
21. Staining Protocol for Adherent Cells

Adherent cells need to be carefully washed to avoid the loss of any cells that round up and come off the plate surface. Loose cells may be harvested from the plate or slide surface and treated as suspension cells, while those remaining adherent to the surface should be washed as adherent cells. If the adherent cells are trypsinized, the loose cells can be recombined with the trypsinized pool; alternatively, the loose cells can be recombined with the adherent portion when the analysis is performed. If growing adherent cells in a plate, the entire plate may be gently spun as part of the wash process to sediment any loose floating cells. Avoid trypsinizing cells prior to labeling with a vital dye, like PI. Cell membranes exposed to trypsin could be transiently permeant to vital dyes for a variable time depending upon the cell line. Cells may be labeled with FLICA before or after trypsinization (Figures 3 and 9).

1. Culture cells in TC-flasks and expose to the experimental conditions.
2. Apoptotic cells may detach and begin to float into the media. Save and spin to pellet and include these cells in the analysis.
3. To stain while adherent, go to Section 22. If suspension cells are needed for analysis, go to Step 4.
4. Trypsinize adherent cells; neutralize with trypsin inhibitor present in cell culture media with 20% FBS; pool cells with any pellets created in Step 2; add 2-5 mL media.
5. Centrifuge at 200 x g for 5 minutes.
6. Remove all but ~100 μ L supernatant. Resuspend cells in 300-500 μ L in cell culture media containing 10-20% FBS. If necessary, count cells and adjust the volume of cell suspension to fit the experiment. Transfer cells into a 15 mL tube.
7. Add 30X FLICA at 1:30.
8. Incubate 30-60 minutes at 37°C, mixing gently every 10 minutes.
9. Wash by adding 2 mL 1X Apoptosis Wash Buffer.
10. Centrifuge at 200 x g for 5 minutes.
11. Aspirate supernatant and resuspend cells in 2 mL 1X Apoptosis Wash Buffer.
12. Incubate 10 minutes at 37°C to allow any unbound FLICA to diffuse out of the cells.
13. Centrifuge at 200 x g for 5 minutes.

Figure 8: Caspase Activity Analyzed with a Fluorescence Plate Reader

Jurkat cells were treated with either DMSO (negative, non-induced cells; left bar) or staurosporine (apoptotic, induced cells; right bar) for 2 hours at 37°C. Cells were labeled with ICT's poly caspase inhibitor reagent, FAM-VAD-FMK (kit #92), for 60 minutes at 37°C. Samples were read on a Molecular Devices Gemini XS 96-well fluorescence plate reader set at 490 nm excitation and 520 nm emission using a 495 cut-off filter. In the induced population, the relative fluorescence units (RFU) of the green fluorescent signal was five times greater than the RFU of the non-induced population (34.11 vs. 6.35). Staurosporine induced poly caspase activity in Jurkat cells. Data courtesy of Dr. Brian Lee, ICT (insert 090309).



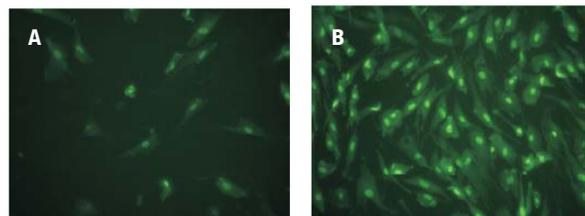
14. Aspirate supernatant and resuspend cells in 1X Apoptosis Wash Buffer. Store cells on ice and protect from light; read within 4 hours. At this point, cells may be analyzed with a fluorescence microscope, flow cytometer, or plate reader.
 - If viewing under a microscope, stain nuclei with Hoechst 33342 according to Step 6 in Section 16. Follow Section 17 for microscopy analysis instructions.
 - If analyzing via flow cytometry, cells do not need to be counted. Follow Section 18 or 19 (Figure 5).
 - If analyzing trypsinized adherent cells (in suspension) with a fluorescence plate reader, follow Section 20.

22. Fluorescence Plate Reader Analysis of Adherent Cells

1. Adherent cells can be grown in a microtiter plate. Use tissue culture plates with clear bottoms and black walls. Culture cells to approximately 90% confluency.
2. Add 30X FLICA at 1:30.
3. Incubate 30-60 minutes at 37°C. Mix gently every 10 minutes to ensure an even distribution of FLICA.
4. Wash by adding ~400 μ L media to each well.
5. Incubate 60 minutes at 37°C to allow any unbound FLICA to diffuse out of the cells.
6. Gently centrifuge the entire plate to sediment any loose floating cells.
7. Aspirate the media. Resuspend with fresh media or PBS.
8. Read plates using a bottom-reading instrument. Set the excitation wavelength at 488 nm and the emission wavelength to 530 nm; use a cut-off filter of 515 nm. FAM-FLICA has an optimal excitation range of 488 - 492 nm and emission range of 515 - 535 nm.

Figure 9: Adherent Corneal Fibroblasts

Normal (A) and keratoconus (B) corneal fibroblasts were treated with 200 μ M H₂O₂ for 1 hour, washed, and allowed to recover for 1-3 hours. The culture media was removed and replaced with ICT's FAM-FLICA caspase-3/7 inhibitor reagent, FAM-DEVD-FMK (kit #94), in cell culture media at 300 μ L/well for 1 hour. The cell layer was washed 3 times with 1X Apoptosis Wash Buffer; 300 μ L was used to keep the cells from drying. Keratoconus corneal fibroblasts treated with H₂O₂ (B) show a significant increase in caspase-3/7 activity compared to normal cells (A). Non-apoptotic cells are dark in background. Data courtesy of Dr. Cristina Kenney, MD, PhD, Dept. of Ophthalmology, University of California, Irvine.



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

23. Staining Protocol for Tissue Sections

See Figure 10.

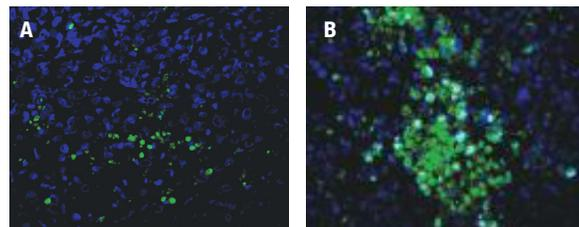
1. Prepare frozen tissues according to the experiment.
2. Dry slides and fix with acetone for 1 minute.
3. Rehydrate slides by washing in TBS-Tween (TBSt) or PBS-Tween (PBSt) for 5 minutes.
4. Wash again in TBSt or PBSt for 5 minutes.
5. Block slides for 20 minutes. Use a blocker such as 20% Aquablock in media with 0.2% Tween.
6. Dilute 150X FLICA stock 1:50 in PBS to form the tissue section staining solution (TSSS). For example, add 50 μ L 150X stock to 2450 μ L PBS (2.5 mL total).
7. Add 50 μ L TSSS per sample (or enough to cover the tissue) and incubate >1 hour protected from light.
8. Wash in TBSt or PBSt for at least 10 minutes.
9. Wash 2 more times in TBSt or PBSt for at least 10 minutes each time.
10. Set slides in a dish containing 1X Apoptosis Wash Buffer.
11. Stain nuclei with Hoechst 33342 (catalog #639) or DAPI (catalog #6244) and apply a coverslip.
12. Store samples at 2-8°C for short-term storage or -20°C for long-term storage (Figure 10).

24. References and Citations

1. Amstad, P. A., et al. Detection of caspase activation *in situ* by fluorochrome-labeled caspase inhibitors. *Biotechniques*. 31, 608-614 (2001).
2. Bedner, E., et al. Activation of caspases measured *in situ* by binding of fluorochrome-labeled inhibitors of caspases (FLICA): correlation with DNA fragmentation. *Exp Cell Res*. 259(1):308-313 (2000).
3. Cascao, R., et al. Caspase-1 is active since the early phase of rheumatoid arthritis. *Ann Rheum Dis*. 70:A2 (2011).
4. Chwa, M., et al. Increased stress-induced generation of reactive oxygen species and apoptosis in human keratoconus fibroblasts. *Invest Ophthalmol Vis Sci*. 47(5): 1902-1910 (2006).
5. Cuevas-Ramos, G., et al. *Escherichia coli* induces DNA damage *in vivo* and triggers genomic instability in mammalian cells. *PNAS*. 107:11537-11542 (2010).
6. Darzynkiewicz, Z., et al. Detection of caspases activation *in situ* by fluorochrome-labeled inhibitors of caspases (FLICA). *Methods Mol Biol*. 203:289-299 (2002).
7. Dunne, A., et al. Inflammasome activation by adenylate cyclase toxin directs Th17 responses and protection against *Bordetella pertussis*. *J Immunol*. 185:1711-1719 (2010).
8. Ekert, P. G., et al. Caspase inhibitors. *Cell Death Differ*. 6:1081-1086 (1999).
9. Evangelisti, C., et al. Preclinical testing of the Akt inhibitor triciribine in T-cell acute lymphoblastic leukemia. *J Cell Physiol*. 226(3):822-31 (2011).
10. Fang, C., et al. The capsid proteins of Aleutian mink disease virus activate caspases and are specifically cleaved during infection. *J Virol*. 84(6):2687-2696 (2010).
11. Fiala, M., et al. HIV-1 induces cardiomyopathy by cardiomyocyte invasion and gp120, Tat and cytokine apoptotic signaling. *Cardiovasc Toxicol*. 4(2):97-107 (2004).
12. Freishtat, R. J., et al. Asthmatic airway epithelium is intrinsically inflammatory and mitotically dyssynchronous. *Am J Respir Cell Mol Biol*. 44:863-869 (2011).
13. Grabarek, J., et al. Use of fluorescently labeled caspase inhibitors as affinity labels to detect activated caspases. *Human Cell*. 15(1):1-12 (2002).
14. Grunewald, S., et al. Caspase activity in human spermatozoa in response to physiological and pathological stimuli. *Fertil Steril* 83:1106-1112 (2005).
15. Kumar, S. Mechanisms mediating caspase activation in cell death. *Cell Death Differ*. 6:1060-1066 (1999).
16. Li, D., et al. Enzymatic dysfunction of mitochondrial complex I of the *Candida albicans* *goal*1 mutant is associated with increased reactive oxidants and cell death. *Eukaryotic Cell*. 10(5):672 (2011).
17. Lopez-Marure, R., et al. Effects of dehydroepiandrosterone on proliferation, migration, and death of breast cancer cells. *Eur J Pharmacol*. 660(2-3):268-74 (2011).
18. Luthra, S. et al. Activation of caspase-8 and caspase-12 pathways by 7-keto-cholesterol in human retinal pigment epithelial cells. *Invest Ophthalmol Vis Sci*. 47:5569-5575 (2006).
19. Mansoor, S. et al. Inhibition of apoptosis in human retinal pigment epithelial cells treated with benzo(e)pyrene, a toxic component of cigarette smoke. *Invest Ophthalmol Vis Sci*. 51:2601-2607 (2010).
20. Olivares-Zavaleta, N., et al. *Chlamydia pneumoniae* inhibits activated human T lymphocyte proliferation by the induction of apoptotic and pyroptotic pathways. *J Immunol*. 186:7120-7126 (2011).
21. Osuka, A., et al. A protective role for inflammasome activation following injury. *Shock*. (2011, Sep 3 Epub ahead of print).
22. Pozarowska, D., et al. Cytometric assessment of cytostatic and cytotoxic effects of topical glaucoma medications on human epithelial corneal line cells. *Cytometry B Clin Cytom*. 78(2):130-7 (2010).
23. Rausch, V., et al. Synergistic activity of sorafenib and sulforaphane abolishes pancreatic cancer stem cell characteristics. *Cancer Res*. 70:5004-5013 (2010).
24. Rotonda, J., et al. The three-dimensional structure of apopain/CPP32, a key mediator of apoptosis. *Nature Struct. Biol*. 3(7):619-625 (1996).
25. Slee, E. A., et al. Serial Killers: ordering caspase activation events in apoptosis. *Cell Death Differ*. 6:1067-1074 (1999).
26. Smoleski, P., et al. Assay of caspase activation *in situ* combined with probing plasma membrane integrity to detect three distinct stages of apoptosis. *J Immunol Methods*. 265(1-2):111-121 (2002).
27. Thomson, L., et al. Cryopreservation-induced human sperm DNA damage is predominantly mediated by oxidative stress rather than apoptosis. *Hum Reprod*. 24:2061-2070 (2009).
28. Thornberry, N. A., et al. 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 (1997).
29. Walker, N. P., et al. Crystal structure of the cysteine protease interleukin-1 β -converting enzyme: a (p20/p10)₂ homodimer. *Cell*. 78:343-352 (1994).
30. Wei, J., et al. Temporal regulation of intracellular organelle homeostasis in T lymphocytes by autophagy. *J Immunol*. 186:5313-5322 (2011).
31. Wilson, K. P., et al. Structure and mechanism of interleukin-1 beta converting enzyme. *Nature*. 370:270-275 (1994).
32. Wolbers, F., et al. Apoptotic cell death kinetics *in vitro* depend on the cell types and the inducers used. *Apoptosis*. 9:385-392 (2004).
33. Yaacob, N. S., et al. Anticancer activity of a sub-fraction of dichloromethane extract of *Strobilanthes crispus* on human breast and prostate cancer cells *in vitro*. *BMC Complement Altern Med*. 10:42 (2010).

Figure 10: Frozen Tumor Sections

A human tumor xenograph was grown subcutaneously in nude mice. Unfixed frozen 10 μ m sections were prepared and stained with ICT's green FLICA poly caspase inhibitor reagent (FAM-VAD-FMK, kit #92) and counterstained with DAPI (blue); total magnification is 400X. A low level of caspase activity was detected in photo A compared with photo B, which reveals extensive green staining in cells throughout the tissue section. Data courtesy of Dr. Rolf Brekken, UTSW.



FOR RESEARCH USE ONLY.

NOT FOR USE IN DIAGNOSTIC PROCEDURES.

Thank you for using FLICA®! If you have any questions, or would like to share your data, please contact us at 1-800-829-3194 or 952-888-8788, or send an email to help@immunochemistry.com.



ImmunoChemistry Technologies, LLC
9401 James Avenue South #155
Bloomington, MN 55431 USA

toll-free: 1-800-829-3194
local: 952-888-8788
fax: 952-888-8988
www.immunochemistry.com