



# Mit-E- $\Psi$ <sup>TM</sup> Mitochondrial Permeability Detection Kit

**Instruction Manual**  
**BML-AK116**

*For research use only*

# MItt-E- $\Psi$ <sup>TM</sup> Mitochondrial Permeability Detection Kit

## BML-AK116

### SECTION 1 Introduction

Detection of the change in mitochondrial permeability provides an early indication of the initiation of cellular apoptosis. This process is typically defined as a collapse in the electrochemical gradient across the mitochondrial membrane, as measured by the change in the membrane potential ( $\Delta\Psi$ ).

Changes in the mitochondrial  $\Delta\Psi$  lead to the insertion of proapoptotic proteins into the membrane and possible oligomerization of BID, BAK, BAX or BAD. This could create pores, which dissipate the transmembrane potential, thereby releasing cytochrome c into the cytoplasm (1-5).

Loss of mitochondrial  $\Delta\Psi$ , indicative of apoptosis, can be detected by a unique fluorescent cationic dye, 5,5',6,6'-tetrachloro-1,1',3,3'-tetraethylbenzamidazolocarboxyanin iodide, commonly known as JC-1 (6). This dye has been incorporated into the user-friendly **MItt-E- $\Psi$ <sup>TM</sup>** kit for the simple and reproducible detection of the change in mitochondrial membrane potential in apoptotic cells.

The structure of Enzo Life Sciences's **MItt-E- $\Psi$ <sup>TM</sup>** reagent allows it to easily penetrate cells and healthy mitochondria. Once inside a healthy non-apoptotic cell, the lipophilic **MItt-E- $\Psi$ <sup>TM</sup>** reagent, bearing a delocalized positive charge, enters the negatively charged mitochondria where it aggregates and fluoresces red (7). These aggregates, first described by Jelley in 1937, are referred to as J-aggregates (8). When the mitochondrial  $\Delta\Psi$  collapses in apoptotic cells, the **MItt-E- $\Psi$ <sup>TM</sup>** reagent no longer accumulates inside the mitochondria. Instead, it is distributed throughout the cell. When dispersed in this manner, the **MItt-E- $\Psi$ <sup>TM</sup>** reagent assumes a monomeric form, which fluoresces green (9). Use of the **MItt-E- $\Psi$ <sup>TM</sup>** kit allows the easy distinction between non-apoptotic red fluorescent cells and apoptotic green fluorescent cells.

The **MItt-E- $\Psi$ <sup>TM</sup>** kit is used in conjunction with your existing apoptosis protocols. Grow your cells in culture and induce apoptosis according to your existing procedure (also reserve a non-induced population of cells as a negative control). Once you have induced apoptosis in your cells, add the 1X **MItt-E- $\Psi$ <sup>TM</sup>** solution to each population and incubate the cells for an additional 15 minutes. During this incubation time, the **MItt-E- $\Psi$ <sup>TM</sup>** reagent will enter each cell and the mitochondria inside. If the cell is not undergoing apoptosis, the mitochondrial  $\Delta\Psi$  will be intact, and the **MItt-E- $\Psi$ <sup>TM</sup>** reagent will aggregate inside the mitochondria and fluoresce red. If the cell

is apoptotic, the mitochondrial  $\Delta\Psi$  will be breaking down, thereby causing the **MItt-E- $\Psi$ <sup>TM</sup>** reagent to be dispersed throughout the entire cell and fluoresce green.

The **MItt-E- $\Psi$ <sup>TM</sup>** reagent excites at 488-490 nm. The monomeric dye structure emits at 527 nm, whereas the J-aggregates in healthy (non-apoptotic) mitochondria emit at 590 nm (7).

The **MItt-E- $\Psi$ <sup>TM</sup>** kit can evaluate apoptosis using three different technologies: flow cytometers; fluorometric plate readers; and fluorescence microscopes.

Following the flow cytometer and fluorescence microscope protocols, each sample to be stained requires only 0.5 mL of 1X **MItt-E- $\Psi$ <sup>TM</sup>** solution (equal to 5  $\mu$ L of 100X **MItt-E- $\Psi$ <sup>TM</sup>** stock). The **MItt-E- $\Psi$ <sup>TM</sup>** Kit will test 100 samples. Following the fluorescence plate reader protocol, each sample requires 1 mL of 1X **MItt-E- $\Psi$ <sup>TM</sup>** solution (equal to 10  $\mu$ L of 100X **MItt-E- $\Psi$ <sup>TM</sup>** stock) and will test 50 samples.

When cells stained with the **MItt-E- $\Psi$ <sup>TM</sup>** reagent are run through a flow cytometer, the instrument will measure apoptosis by monitoring the amount of red fluorescence in each region. Healthy cells, which fluoresce red, will appear in R2. As the mitochondrial  $\Delta\Psi$  collapses and cells enter apoptosis, the amount of red fluorescence will drop. An increasing number of cells will fall into R3 corresponding to a loss of red fluorescence as the dispersed **MItt-E- $\Psi$ <sup>TM</sup>** dye converts to a monomeric form and fluoresces green (see Section 16 for sample data).

When cells stained with the **MItt-E- $\Psi$ <sup>TM</sup>** reagent are analyzed with a fluorescence plate reader, the instrument will measure apoptosis by monitoring the amount of red fluorescence. Healthy cells will give a high level of red fluorescence emission. As the mitochondrial  $\Delta\Psi$  collapses, indicating apoptosis, an increasing number of cells will emit less red fluorescence as the dispersed **MItt-E- $\Psi$ <sup>TM</sup>** dye converts to a green monomeric form (see Section 19 for sample data).

Looking at the cells under a fluorescence microscope, non-apoptotic cells will appear to have red fluorescent spots (the **MItt-E- $\Psi$ <sup>TM</sup>** dye aggregates) within healthy mitochondria. In contrast, apoptotic cells will appear mostly green. Cells still undergoing apoptosis will contain fewer and fewer red dye aggregate spots in the mitochondria, and the entire cell will appear more

and more green as the mitochondrial  $\Delta\Psi$  disintegrates, thereby dispersing the MIt-E- $\Psi^{\text{TM}}$  dye

## SECTION 2 References

1. S. Desagher, *et al. J. Cell Biol.* 1999 **144**: 891.
2. M.Narita, *et al. Proc. Natl. Acad. Sci. USA* 1998 **95**:14681.
3. G.Basanez, *et al. Proc. Natl. Acad. Sci. USA* 1999 **96**:5492.
4. X.Luo, *et al. Cell* 1998 **94**:481.
5. B.Antonsson, *et al. Biochem J.* 2000 **345**:271.
6. S.T.Smiley, *et al. Proc. Natl. Acad. Sci. USA* 1991 **88**:3671.
7. A.Cossarizza, *et al. Biochem. Biophys. Res. Commun.* 1993 **197**: 40.
8. E.Jelley. *Nature* 1939 **139**:631.
9. M.Reers, *et al. Biochemistry* 1991 **30**: 4480.

**PLEASE READ ENTIRE BOOKLET BEFORE PROCEEDING WITH THE ASSAY. CAREFULLY NOTE THE HANDLING AND STORAGE CONDITIONS OF EACH KIT COMPONENT. PLEASE CONTACT TECHNICAL SUPPORT FOR ASSISTANCE IF NECESSARY.**

## SECTION 3 Components

**4x KI-162** MIt-E- $\Psi^{\text{TM}}$  Reagent, lyophilized  
**1x KI-163** 10X Assay Buffer (60 mL)

Assay Manual with protocols for 3 applications: Flow Cytometer; Fluorometer; and Fluorescence Microscope

### Storage and Shelf-Life

- Store the unopened kit (and each unopened component) at 2°C to 8°C until the expiration date.
- Once reconstituted, store the 100X MIt-E- $\Psi^{\text{TM}}$  stock at -20°C for 6 months. It may be frozen twice.

Once diluted, store the 1X assay buffer at 2°C to 8°C up to 7 days.

### Safety Information

- Use gloves while handling the MIt-E- $\Psi^{\text{TM}}$  reagent, and the 10X assay buffer.
- Dispose of all liquid components down the sink and flush with copious amounts of water. Solid components may be disposed of in regular trash.

## SECTION 4 Other Materials

**Depending on experimental protocol used, the following materials may be needed.**

- Pipette(s) capable of dispensing at 10  $\mu\text{L}$ , 500  $\mu\text{L}$ , 1 mL
  - Dimethyl Sulfoxide (DMSO) 125 or 500  $\mu\text{L}$
  - Vortexer
  - Amber vials or polypropylene tubes for storage at -20°C
  - 37°C CO<sub>2</sub> incubator
  - Flow cytometer with excitation between 488 - 490nm, and emission at 527 and 590 nm.
  - 96-well fluorescence plate reader with excitation between 488 -490 nm, and emission at 590 nm (dual emission at 527 nm and 590 - 600 nm is best), with endpoint reading and black round or flat bottom 96-well microtiter plates.
  - Fluorescence microscope with broad band path filters and slides.
  - 96-well microtiter plates, black
- Cultured cells and media
  - Protocol and reagents to induce apoptosis
  - 15 mL polystyrene centrifuge tube (1 per sample)
  - Microcentrifuge at 13,000 X g
  - Clinical centrifuge at 400 – 1,000 X g.

## SECTION 5 Overview of the Mīt-E-Ψ™ Protocol

Staining cells with Mīt-E-Ψ™ takes only about 15 minutes. However, the Mīt-E-Ψ™ kit is used with living cells, which may take several hours to prepare. In addition, once the cells are grown, they must be induced to undergo apoptosis, which may also take several hours to complete. As the 1X Mīt-E-Ψ™ solution must be used immediately, prepare the Mīt-E-Ψ™ reagents at the end of your apoptosis induction process.

**Culture cells to a density optimal for apoptosis induction according to your specific induction protocol, but not to exceed 10<sup>6</sup> cells/mL.**

1. At the same time, culture a non-induced negative control cell population (at the same density as the induced population).
2. Induce apoptosis following your protocol (2 sample protocols are mentioned in Section 6).
3. Prepare 1X assay buffer and warm to 37°C (Section 7).
4. Prepare 100X Mīt-E-Ψ™ stock (Section 8).
5. Prepare 1X Mīt-E-Ψ™ solution (Section 9 or 11).
6. Stain cells with 1X Mīt-E-Ψ™ solution (Section 12, 17, 20, or 22).
7. Analyze data (Section 13, 18, or 22).

## SECTION 6 Induction of Apoptosis

The Mīt-E-Ψ™ kit works with your current apoptosis protocols. 2 examples of protocols to induce apoptosis in suspension culture are: a) treating Jurkat cells with 2 μg/mL camptothecin (or 1 μM staurosporine) for 3 hours; and b) treating HL-60 cells with 4 μg/mL camptothecin (or 1 μM staurosporine) for 4 hours.

## SECTION 7 Preparation of 1X Assay Buffer

The assay buffer is formulated for use as reaction buffer, and for washing the cells. It is supplied as a 10X concentrate which must be diluted to 1X with DI H<sub>2</sub>O prior to use.

1. If necessary, gently warm the 10X concentrate to completely dissolve any salt crystals that may have come out of solution.
2. Add the entire bottle (60 mL) of 10X assay buffer to 540 mL of DI H<sub>2</sub>O. **Or**, if not using the entire bottle, dilute the 10X assay buffer 1:10 in DI H<sub>2</sub>O. For example, add 10 mL 10X assay buffer to 90 mL DI H<sub>2</sub>O.
3. Stir the solution for at least 5 minutes.
4. If using the 1X assay buffer the same day it was prepared, warm it to 37°C prior to

use (it is used in Sections 9, 11, 12, 17, 20, and 21). **Or**, if not using the 1X assay buffer the same day it was prepared, store it at 2°C to 8°C up to 7 days.

## SECTION 8 Reconstitution of the 100X Mīt-E-Ψ™ Stock

The Mīt-E-Ψ™ dye reagent is supplied as a highly concentrated lyophilized powder. It must first be reconstituted, forming a 100X stock concentrate, and then diluted 1:100 to form a final 1X working solution. The 1X working solution must be prepared immediately prior to use; however, the reconstituted 100X stock can be stored at –20°C for 6 months, and used twice during that time.

**The newly reconstituted 100X Mīt-E-Ψ™ stock must be used or frozen immediately after it is prepared and protected from light during handling.**

1. Reconstitute each vial of the Mīt-E-Ψ™ reagent with 125 μL DMSO at room temperature (RT), forming a 100X stock.
2. Re-cap each vial and invert it several times to fully dissolve the Mīt-E-Ψ™ dye reagent.
3. Immediately use the 100X stock by diluting it to 1X (Section 9).
4. **Or**, aliquot and store it at –20°C (Section 10).

## SECTION 9 Preparation of 1X Mīt-E-Ψ™ Solution for Immediate Use

Using the freshly reconstituted 100X Mīt-E-Ψ™ stock, prepare the 1X working strength Mīt-E-Ψ™ solution by diluting the stock 1:100 with 37°C 1X assay buffer (or, you may substitute your own cell culture media, warmed to 37°C, in place of the 1X assay buffer). Each sample to be stained requires only 0.5 mL of 1X Mīt-E-Ψ™ solution (equal to 5 μL of 100X Mīt-E-Ψ™ stock).

**The 1X working strength Mīt-E-Ψ™ solution must be used immediately – prepare it as apoptosis induction is completed.**

1. Warm the 1X assay buffer (or cell culture media) to 37°C.
2. Add 125 μL of the 100X Mīt-E-Ψ™ stock from each vial to 12.375 mL of the 37°C 1X assay buffer or cell culture media. **Or**, add 500 μL (the contents of all 4 vials) of the 100X Mīt-E-Ψ™ stock to 49.5 mL of the 37°C 1X assay buffer or cell culture media. **Or**, if not using the entire vial of 100X Mīt-E-Ψ™ stock, dilute it 1:100 in 37°C 1X assay buffer or cell culture media. For example, add 10 μL of 100X Mīt-

- E-Ψ™ stock to 990 μL of 1X assay buffer or cell culture media.
3. Vortex the 1X working strength Mīt-E-Ψ™ solution thoroughly.
  4. If particulate matter is present, the solution should be clarified by centrifugation at 13,000 X g in a microcentrifuge for 3 minutes, or 15 minutes in a clinical centrifuge at 1,000 X g at RT. This is especially important when using a fluorometer.

**Clarification of the 1X solution is absolutely necessary when performing the fluorometer protocol. Any free dye aggregates in the solution will interfere with the OD reading, leading to a falsely increased reading of red fluorescence. (The aggregates do not interfere with a flow cytometer, and are easily identifiable under a fluorescence microscope.)**

5. If centrifuged, transfer the clarified supernatant to a clean tube and discard particulates.
6. Go on to the staining protocol (Section 12, 17, 20, or 21).

**To avoid photo-bleaching and degradation of the dye, protect the Mīt-E-Ψ™ dye reagent from light while handling.**

## **SECTION 10 Storage of 100X Mīt-E-Ψ™ Stock for Future Use**

If not all of the 100X Mīt-E-Ψ™ stock will be used the same time it is reconstituted, the unused portion may be stored at -20°C for 6 months. During that time, the 100X Mīt-E-Ψ™ stock may be thawed and used twice. After the second thaw, discard any remaining 100X Mīt-E-Ψ™ stock. If you anticipate using it more than twice, make small aliquots and store in amber vials or polypropylene tubes.

Following the flow cytometer and fluorescence microscope protocols, each sample to be stained requires only 0.5 mL of 1X Mīt-E-Ψ™ solution (equal to 5 μL of 100X Mīt-E-Ψ™ stock). Following the fluorescence plate reader protocol, each sample requires 1 mL of 1X Mīt-E-Ψ™ solution (10 μL of 100X Mīt-E-Ψ™ stock). When needed, prepare frozen Mīt-E-Ψ™ solution according to directions in Section 11.

## **SECTION 11 Preparation of 1X Mīt-E-Ψ™ Solution from a Frozen Aliquot**

If some of the 100X Mīt-E-Ψ™ stock was previously reconstituted and then stored at -20°C, it may be used 2 more times within 6 months.

1. Thaw the 100X Mīt-E-Ψ™ stock and protect from light.

2. Once the aliquot has become liquid, dilute the 100X Mīt-E-Ψ™ stock 1:100 in 37°C 1X assay buffer or cell culture media. For example, mix 10 μL of 100X Mīt-E-Ψ™ stock with 990 μL of 37°C 1X assay buffer or cell culture media.
3. If the 100X Mīt-E-Ψ™ stock was frozen immediately after reconstitution and was never thawed, return it to the freezer. If the stock has been thawed once before, discard it.
4. Vortex the 1X working strength Mīt-E-Ψ™ solution thoroughly.
5. If particulate matter is present, the solution must be clarified by centrifugation at 13,000 X g in a microcentrifuge for 3 minutes, or 15 minutes in a clinical centrifuge at 1,000 X g at RT.

**Clarification of the 1X solution is especially important when performing the fluorometer protocol. Any free dye aggregates in the solution will interfere with the OD reading, leading to a falsely increased reading of red fluorescence. The aggregates do not interfere with a flow cytometer, and are easily identifiable under a fluorescence microscope.)**

6. If centrifuged, transfer the clarified supernatant to a clean tube and discard particulates.
7. Go on to the staining protocol (Section 12, 17, 20, or 21).

## SECTION 12 Flow Cytometry Staining Protocol

Following the flow cytometer protocols, each sample to be stained requires only 0.5 mL of 1X MIt-E- $\Psi^{\text{TM}}$  solution (equal to 5  $\mu\text{L}$  of 100X MIt-E- $\Psi^{\text{TM}}$  stock). The MIt-E- $\Psi^{\text{TM}}$  Kit will test 100 samples.

1. As discussed in Section 6, culture cells to a density optimal for apoptosis induction according to your specific induction protocol.
2. Induce apoptosis following your protocol.
3. At the same time, culture a non-induced negative control cell population.
4. After the induction process, transfer 0.5 mL of each cell suspension into a sterile 15 mL polystyrene centrifuge tube.

**Cell density in the cell culture flasks should not exceed  $10^6$  cells per mL. Cells cultivated in excess of this concentration may begin to naturally enter apoptosis. Optimal cell concentration will vary depending on the cell line used. Density can be determined by counting cell populations on a hemocytometer.**

5. Centrifuge cells at 400 X g for 5 minutes at RT.
6. Carefully remove and discard the supernatant.
7. Gently vortex the cell pellets, or use a pipette, to disrupt any cell-to-cell clumping.
8. Resuspend cells in 0.5 mL of RT 1X MIt-E- $\Psi^{\text{TM}}$  solution.
9. Incubate the cells (which are now being stained with the MIt-E- $\Psi^{\text{TM}}$  dye reagent) at 37°C for 10-15 minutes in a CO<sub>2</sub> incubator.
10. Warm the 1X assay buffer to 37°C (Section 7).
11. Add 2 mL of 1X assay buffer to each tube.
12. Mix each tube.
13. Centrifuge the stained cells at 400 X g for 5 minutes at RT.
14. Carefully remove and discard supernatant.
15. Gently vortex the pellet to disrupt any cell-to-cell clumping.
16. Resuspend the cells in 1 mL of 1X assay buffer.
17. Centrifuge the stained cells at 400 X g for 5 minutes at RT.
18. Carefully remove and discard supernatant.
19. Gently vortex the pellet to disrupt any cell-to-cell clumping.
20. Resuspend the cells in 1 mL of 1X assay buffer.
21. Centrifuge the stained cells at 400 X g for 5 minutes at RT.
22. Carefully remove and discard supernatant.
23. Gently vortex the pellet to disrupt any cell-to-cell clumping.

24. Resuspend the cell pellet in 0.5 mL of 1X assay buffer.
25. Analyze cells by flow cytometry (Sections 13-16).

## SECTION 13 Flow Cytometer Set Up

Healthy cells containing the aggregated MIt-E- $\Psi^{\text{TM}}$  dye reagent within their mitochondria (which fluoresces red) can be detected in the FL2 channel. Apoptotic cells, which contain non-aggregated green MIt-E- $\Psi^{\text{TM}}$  monomers, can be detected in the FL1 channel.

## SECTION 14 Single-Parameter Analysis Using a Flow Cytometer

1. Generate a log FL2 (X-axis) versus cell count.
2. Adjust the FL2 PMT voltage to allow the peak to fall within the third log decade when running the negative control sample.
3. When running the induced positive samples, use the same adjusted PMT voltage as was determined for the negative control.
4. Observe the mean fluorescence of both the apoptotic positive cell population, as well as the mean fluorescence of the treated negative control cell population.

**The apoptotic cell population presents a lower red fluorescence signal intensity (FL2 axis) than the negative control population.**

## SECTION 15 Multi-Parameter Analysis Using a Flow Cytometer

1. Create a log FL1 (X-axis) versus log FL2 (Y-axis) scatter plot.
2. Add two regions, R2 and R3, as shown in Figure 1.
3. Run the negative control sample first (non-induced cells). Adjust the FL1 and FL2 PMT voltages so that the majority of the cell population falls within the upper right hand region. The peak of the dual fluorescent population should fall within the second and third log decade scale of FL1 (X-axis) and FL2 (Y-axis) as seen in Figure 1.
4. Adjust R2 so that greater than 95% of the dual fluorescent cell population falls within this region. The number of cells falling into this region will vary depending on the condition of the culture and cell type.
5. Next, adjust R3 so that it falls directly below R2, as in Figure 1.
6. Using the same PMT settings established for the non-induced negative control

sample in Step 3, run the induced positive sample. If a change in mitochondrial  $\Delta\Psi$  has occurred, an increase in the number of cells falling in R3 is observed as shown in Figure 2. **This reflects a reduction in red fluorescence.**

7. If the induced sample exhibits only a minimal change in red emission, increase the FL2-FL1 compensation and repeat Steps 3-6.

## SECTION 16 Flow Cytometry Sample Data

When cells stained with MIt-E- $\Psi^{\text{TM}}$  are run through a flow cytometer, the instrument will measure apoptosis by monitoring the amount of red fluorescence in each region. Healthy cells, which fluoresce red, will appear in R2. As the mitochondrial  $\Delta\Psi$  collapses and cells enter apoptosis, the amount of red fluorescence will drop. An increasing number of cells will fall into R3 corresponding to a loss of red fluorescence as the dispersed MIt-E- $\Psi^{\text{TM}}$  dye converts to a monomeric form and fluoresces green.

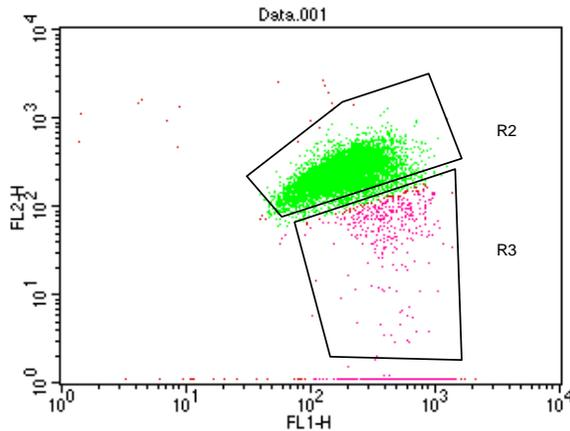


Figure 1: Negative cell control, Log FL1 (X-axis) and log FL2 (Y-axis) R2 = 91.07% R3 = 8.08%.

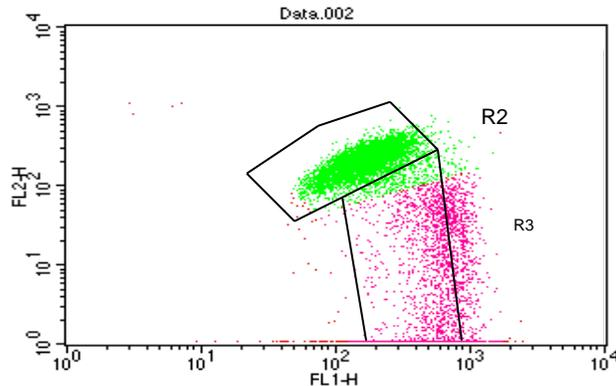


Figure 2: Positive cell control, Log FL1 (X-axis) and log FL2 (Y-axis) R2 = 60.93% R3 = 38.13%.

In Figures 1 and 2, cells were analyzed in a FACS Caliber Becton Dickinson Flow Cytometer. Jurkat cells were either treated with DMSO (negative, non-induced cells) or with staurosporine (apoptotic, induced cells) for 3 hours at 37°C and then labeled with MIt-E- $\Psi^{\text{TM}}$  for 15 minutes. Collapse of the mitochondrial  $\Delta\Psi$  is indicated by an increase in the number of cells falling into R3 corresponding to a loss of red fluorescence, indicative of the onset of apoptosis. (In these figures of flow cytometer data, non-apoptotic cells are in R2, and apoptotic cells are in R3.)

## SECTION 17

### 96-Well Fluorescence Spectroscopy Staining Protocol

Following the fluorescence plate reader protocol, each sample requires 1 mL of 1X MIt-E- $\Psi^{\text{TM}}$  solution (equal to 10  $\mu\text{L}$  of 100X MIt-E- $\Psi^{\text{TM}}$  stock). The MIt-E- $\Psi^{\text{TM}}$  Kit will test 50 samples.

1. As discussed in Section 6, culture cells to a density optimal for apoptosis induction according to your specific induction protocol.

**Cell density in the cell culture flasks should not exceed  $10^6$  cells per mL. Cells cultivated in excess of this concentration may begin to naturally enter apoptosis. Optimal cell concentration will vary depending on the cell line used. Density can be determined by counting cell populations on a hemocytometer.**

2. Induce apoptosis following your protocol.
3. At the same time, culture an equal volume of non-induced cells for a negative control cell population. Make sure that both tubes of cells contain similar quantities of cells.
4. Pellet cells by centrifugation at 400 X g for 5 minutes at RT.

**When concentrated, enough cells should have been grown to form a 0.5 mL concentrated pool between 1 - 2 X  $10^6$  cells/mL.**

5. Carefully remove and discard the supernatants.
6. Resuspend at between 0.5 – 2 X  $10^6$  cells in 1 mL of 1X working strength MIt-E- $\Psi^{\text{TM}}$  solution (Sections 9 and 11).
7. Gently vortex the cell pellets, or use a pipette, to disrupt any cell-to-cell clumping.
8. Incubate the cells (which are now being stained with the MIt-E- $\Psi^{\text{TM}}$  dye reagent) at 37°C for 10-15 minutes in a CO<sub>2</sub> incubator.
9. Warm the working strength 1X assay buffer to 37°C (Section 7).
10. Add 2 mL of 1X assay buffer to each tube.
11. Mix each tube.
12. Centrifuge the stained cells at 400 X g for 5 minutes at RT.
13. Carefully remove and discard supernatants.
14. Gently vortex the pellets to disrupt any cell-to-cell clumping.
15. Resuspend the cells in 1 mL of 1X assay buffer.
16. Take out a small aliquot to determine the concentration of both the induced and non-induced cell populations.
  - a. Remove a 50  $\mu\text{L}$  aliquot of each cell population.
  - b. Add to 450  $\mu\text{L}$  PBS (forming a 1:10 dilution of each).

- c. Count the cells (a hemocytometer may be used).
- d. After counting, compare the density of each. The non-induced population may have more cells than the induced population, as some induced cells may be lost during the apoptotic process. If necessary, adjust the volume of the induced cell suspension to match that of the non-induced suspension (Step 20).

17. Centrifuge the remaining stained cells at 400 X g for 5 minutes at RT.
18. Carefully remove and discard supernatants.
19. Gently vortex the pellets to disrupt any cell-to-cell clumping.
20. Adjust the volume of the induced cell suspension to match that of the non-induced suspension.

**At least 0.5 X  $10^6$  cells/mL are required to generate an adequate signal using a fluorometer.**

- a. Resuspend the non-induced cell pellets in 500  $\mu\text{L}$  - 1 mL of 1X assay buffer to produce a cell suspension at least 5 X  $10^5$  cells/mL. The volume may vary depending upon cell density.
  - b. Resuspend the induced cell population in a volume of 1X assay buffer to yield the same concentration of cells as the non-induced cell suspension.
21. For each sample to be tested, dispense 100  $\mu\text{L}$  into each of 2 wells in a black round or flat bottom 96-well microtiter plate.

**At least 5 X  $10^4$  cells/well are required to generate an adequate signal using a fluorometer, with best results obtained using 1 X  $10^5$  cells/well. The 96-well microtiter plate used to analyze the cells must be black. A clear plate will interfere with the OD readings.**

22. Analyze cells using a fluorescence plate reader (Section 18).

## SECTION 18

### 96-Well Fluorescence Plate Reader Set Up

1. Set the plate reader to perform an endpoint read.
2. Set the excitation wavelength at 488 - 490 nm
3. Set the emission wavelengths to 527 nm for green fluorescence and 590 - 600 nm for red fluorescence. If your plate reader cannot read dual emission wavelengths at the same time, use the red fluorescence setting of 590 - 600 nm to perform the analysis.
4. Read the samples.

## SECTION 19 96-Well Fluorescence Spectroscopy Sample Data

When cells stained with MIt-E- $\Psi^{\text{TM}}$  are analyzed with a fluorescence plate reader, the instrument will measure the amount of red fluorescence. Healthy cells will emit a high level of red fluorescence; apoptotic cells will generate a lower level of red fluorescence.

By comparing the average 590-600 nm emission levels in stimulated versus non-stimulated sample wells, loss of mitochondrial  $\Delta\Psi$  can be monitored. As the mitochondrial  $\Delta\Psi$  collapses (indicating apoptosis) and the dispersed MIt-E- $\Psi^{\text{TM}}$  dye converts to a green monomeric form, more and more cells will lose red fluorescence (Figures 3 and 4).

Using the dual fluorescence characteristic of the dye, the changes in the mitochondrial  $\Delta\Psi$  can be most accurately assessed by comparing the ratio of 590-600 nm (red) / 527 nm (green) emissions. When apoptosis is induced, the red/green emission ratio drops compared to the negative (non-stimulated) control wells (the red emission level decreases and the green emission level increases). This drop corresponds to a reduction in the number of healthy mitochondria able to maintain the negative potential necessary to concentrate the MIt-E- $\Psi^{\text{TM}}$  dye in the red aggregate form.

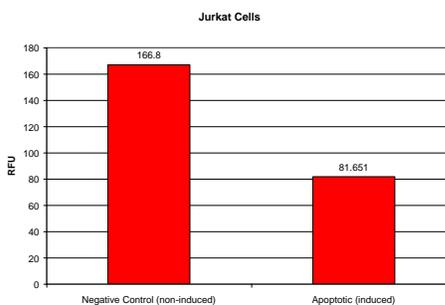


Figure 3: Red fluorescence of Jurkat cells

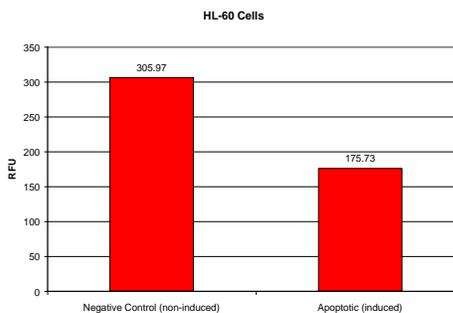


Figure 4: Red fluorescence of HL-60 cells

In Figures 3 and 4, the cells were either treated with DMSO (negative, non-induced cells) or with staurosporine (apoptotic, induced cells) for 4 hours and then labeled with the 1X MIt-E- $\Psi^{\text{TM}}$  solution for 15 minutes. The samples were then read on a 96-well fluorescence plate reader using the settings described above. As the mitochondrial  $\Delta\Psi$  collapses, indicating apoptosis, the amount of red fluorescence drops by 51% in the Jurkat cells and 43% in HL-60 cells.

**If there are free aggregates of the MIt-E- $\Psi^{\text{TM}}$  dye in solution with the cells, they will fluoresce red and interfere with the data, leading to a falsely increased reading of red fluorescence (Sections 9 and 11).**

## **SECTION 20**

### **Fluorescence Microscopy Staining Protocol for Adherent Cells**

Following the fluorescence microscopy protocol, each sample to be stained requires only 0.5 mL of 1X MIt-E- $\Psi$ <sup>TM</sup> solution (equal to 5  $\mu$ L of 100X MIt-E- $\Psi$ <sup>TM</sup> stock). The MIt-E- $\Psi$ <sup>TM</sup> Kit will test 100 samples.

1. Culture cells on a sterile coverslip or chamber slide to a cell density optimal for apoptosis induction according to your specific induction protocol.

**Cell density should not exceed the threshold where cell sloughing occurs.**

2. As discussed in Section 6, induce apoptosis following your protocol.
3. At the same time, culture an equal volume of non-induced cells for a negative control cell population.
4. Make sure that both the negative control and induced positive cell populations contain similar quantities of cells.
5. Remove the media from induced and non-induced monolayer cultures.
6. Add enough 1X MIt-E- $\Psi$ <sup>TM</sup> solution (Sections 9 and 11) to cover the cells on the slide.
7. Incubate the cells (which are now being stained with the MIt-E- $\Psi$ <sup>TM</sup> dye reagent) at 37°C for 15 minutes in a CO<sub>2</sub> incubator.
8. Warm the 1X assay buffer to 37°C (Section 7).
9. Carefully remove and discard staining media.
10. Wash the monolayer cultures with 1 to 2 mL of 1X assay buffer.
11. Discard wash.
12. Add a drop of 1X assay buffer plus coverslip.
13. Examine using a fluorescence microscope (Section 22).

## **SECTION 21**

### **Fluorescence Microscopy Staining Protocol for Suspension Cells**

Following the fluorescence microscope protocol, each sample to be stained requires only 0.5 mL of 1X MIt-E- $\Psi$ <sup>TM</sup> solution (equal to 5  $\mu$ L of 100X MIt-E- $\Psi$ <sup>TM</sup> stock). The MIt-E- $\Psi$ <sup>TM</sup> Kit will test 100 samples.

1. Culture cells to a cell density optimal for apoptosis induction according to your specific induction protocol.

**Optimal cell densities will vary with the cell line; cell concentrations may be determined using a hemo-cytometer.**

2. As discussed in Section 6, induce apoptosis following your protocol.

3. At the same time, culture an equal volume of non-induced cells for a negative control cell population.
4. Make sure that both the negative control and induced positive cell populations contain similar quantities of cells.
5. Once induced, count cells (a hemocytometer may be used).
6. Transfer 0.5 X 10<sup>6</sup> cells to a centrifuge tube.
7. Centrifuge cells at 400 X g for 5 minutes at RT.
8. Carefully remove and discard the supernatants.
9. Gently vortex the cell pellets, or use a pipette, to disrupt any cell-to-cell clumping.
10. Resuspend cells in 0.5 mL 1X MIt-E- $\Psi$ <sup>TM</sup> solution (Sections 9 and 11).
11. Incubate the cells (which are now being stained with the MIt-E- $\Psi$ <sup>TM</sup> dye reagent) at 37°C for 10 to 15 minutes in a CO<sub>2</sub> incubator.
12. Warm the 1X assay buffer to 37°C (Section 7).
13. Resuspend cells in 2 mL 1X assay buffer.
14. Centrifuge cells at 400 X g for 5 minutes at RT.
15. Carefully remove and discard the supernatants.
16. Gently vortex the cell pellets to disrupt any cell-to-cell clumping.
17. Resuspend cells in 1 mL of 1X assay buffer.
18. Centrifuge the stained cells at 400 X g for 5 minutes at RT.
19. Carefully remove and discard the supernatants.
20. Gently vortex the cell pellets to disrupt any cell-to-cell clumping.
21. Resuspend cells in 1 mL of 1X assay buffer.
22. Centrifuge the stained cells at 400 X g for 5 minutes at RT.
23. Carefully remove and discard the supernatants.
24. Resuspend cells in 0.5 mL 1X assay buffer.
25. Examine a drop of cell suspension under a cover slip using a fluorescence microscope (Section 22).

## **SECTION 22**

### **Fluorescence Microscope Set Up**

This scope should contain a long band path filter (Ex 490 nm, Em >510 nm) capable of detecting both fluorescein and rhodamine fluorescence.

## **DISCLAIMER**

**NOTE: THE APPLICATION EXAMPLES, DESCRIBED ABOVE, ARE INTENDED ONLY AS GUIDELINES. THE OPTIMAL CONCENTRATIONS OF SUBSTRATES AND INHIBITORS, ASSAY VOLUMES, BUFFER COMPOSITION, AND OTHER EXPERIMENTAL CONDITIONS MUST BE DETERMINED BY THE INDIVIDUAL USER. NO WARRANTY OR GUARANTEE OF PARTICULAR RESULTS, THROUGH THE USE OF THESE PROCEDURES, IS MADE OR IMPLIED.**

## **USE OF PRODUCT**

This product contains research chemicals. As such, they should be used and handled only by or under the supervision of technically qualified individuals. This product is not intended for diagnostic or human use.

## **WARRANTY**

Enzo Life Sciences International, Inc makes no warranty of any kind, expressed or implied, which extends beyond the description of the product in this brochure, except that the material will meet our specifications at the time of delivery. Enzo Life Sciences International, Inc makes no guarantee of results and assumes no liability for injuries, damages or penalties resulting from product use, since the conditions of handling and use are beyond our control.

**www.enzolifesciences.com**  
*Enabling Discovery in Life Science®*

---

**Global Headquarters**

**Enzo Life Sciences Inc.**

10 Executive Blvd

Farmingdale, NY 11735

(p) 1-800-942-0430

(f) 1-631-694-7501

(e) [info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**Enzo Life Sciences (ELS) AG**

Industriestrasse 17, Postfach

CH-4415 Lause / Switzerland

(p) +41/0 61 926 89 89

(f) +41/0 61 926 89 79

(e) [info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

Please visit our website at [www.enzolifesciences.com](http://www.enzolifesciences.com) for additional contact information.