

ab65310

**Mitochondrial
Transmembrane
Potential Apoptosis
Detection Kit**

Instructions for Use

For the rapid, sensitive and accurate measurement of apoptosis in various samples.

This product is for research use only and is not intended for diagnostic use.

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

Table of Contents

1.	Overview	3
2.	Protocol Summary	4
3.	Components and Storage	5
4.	Assay Protocol	6
5.	Data Analysis	8
6.	Troubleshooting	9

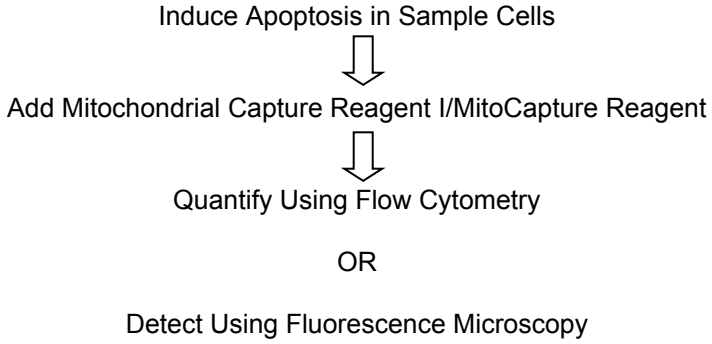
1. Overview

Disruption of the mitochondrial transmembrane potential is one of the earliest intracellular events that occur following induction of apoptosis. Abcam's Mitochondrial Transmembrane Potential Apoptosis Detection Kit provides a simple, fluorescent-based method for distinguishing between healthy and apoptotic cells by detecting the changes in the mitochondrial transmembrane potential.

The kit utilizes Mitochondrial Capture Reagent I/MitoCapture™, a cationic dye that fluoresces differently in healthy vs apoptotic cells. In healthy cells, Mitochondrial Capture Reagent I/MitoCapture accumulates and aggregates in the mitochondria, giving off a bright red fluorescence. In apoptotic cells, Mitochondrial Capture Reagent I/MitoCapture cannot aggregate in the mitochondria due to the altered mitochondrial transmembrane potential, and thus it remains in the cytoplasm in its monomer form, fluorescing green.

The fluorescent signals can be easily detected by fluorescence microscopy using a band-pass filter (detects FITC and rhodamine) or analyzed by flow cytometry using FITC channel for green monomers (Ex/Em = 488/530 ± 30 nm) and (optional) PI channel for red aggregates (Em = 488/590 ± 42 nm).

2. Protocol Summary



3. Components and Storage

A. Kit Components

Item	Quantity (25 assays)	Quantity (100 assays)
Mitochondrial Capture Reagent I/MitoCapture Reagent	25 μ L	100 μ L
Binding Buffer III/Incubation Buffer	50 mL	2 x 100 mL

Store MitoCapture at -20°C. Avoid freeze-thaw cycles. Protect from light. Store Binding Buffer III/Incubation Buffer at +4°C after opening. All reagents are stable for 1 year.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Flow Cytometer or Fluorescence Microscope
- Glass slides and coverslips

4. Assay Protocol

1. Incubation of cells with Mitochondrial Capture Reagent I/MitoCapture Reagent:

- a) Induce apoptosis by desired methods. Concurrently incubate a control culture *without* induction.
- b) Collect $\sim 1 \times 10^6$ cells per sample by centrifugation at 500 x g for 5 minutes.
- c) Aliquot enough Binding Buffer III/Incubation Buffer for the number of assays to be performed (total 2 ml for each assay) and pre-warm to 37°C before use. Dilute Mitochondrial Capture Reagent I/MitoCapture Reagent immediately prior to use: Dilute 1 μ l Mitochondrial Capture Reagent I/MitoCapture to 1 ml in pre-warmed Binding Buffer III/Incubation Buffer for each assay. Vortex the solution.

Note:

Mitochondrial Capture Reagent I/MitoCapture is poorly soluble in aqueous solutions. To remove particles (optional), centrifuge the dye solution for 1 minute at 13,000 x g and carefully transfer the supernatant without disturbing pelleted debris.

- d) Re-suspend cells in 1 ml of the diluted Mitochondrial Capture Reagent I/MitoCapture solution.

- e) Incubate at 37°C in a 5% CO₂ incubator for 15-20 min.
- f) Centrifuge cells at 500 x g and discard supernatant.
- g) Re-suspend in 1 ml of the pre-warmed Binding Buffer III/Incubation Buffer.

Proceed to Step 2 or 3 below depending on method of analysis.

2. Quantification by Flow Cytometry:

Analyze cells immediately following step 1.g by flow cytometry. Mitochondrial Capture Reagent I/MitoCapture monomers in apoptotic cells are detectable in the FITC channel (usually FL1) showing diffused green fluorescence. Mitochondrial Capture Reagent I/MitoCapture aggregates in healthy cells are detectable in the PI channel (usually FL2) showing punctate red fluorescence.

3. Detection by Fluorescence Microscopy:

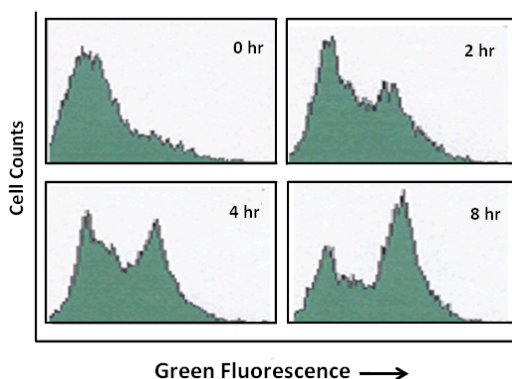
- a) Place the cell suspension from Step 1.g on a glass slide. Cover the cells with a glass coverslip.

For analyzing **adherent cells**, grow cells on a coverslip and perform the entire procedure directly on the coverslip in culture dish. Following incubation (1.e), invert coverslip on a glass slide.

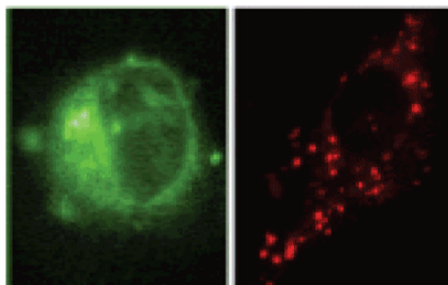
- b) Observe cells immediately under a fluorescence microscope using a band-pass filter (detects fluorescein and rhodamine).

Mitochondrial Capture Reagent I/MitoCapture that has aggregated in the mitochondria of healthy cells fluoresces red. In apoptotic cells, Mitochondrial Capture Reagent I/MitoCapture cannot accumulate in mitochondria, it remains as monomers in the cytoplasm, and fluoresces green.

5. Data Analysis



Apoptosis was induced in Jurkat cells by 2 μ M camptothecin for various hours as indicated. Cells were incubated with Mitochondrial Capture Reagent I/MitoCapture for 15 minutes. Results were analyzed by Flow Cytometry using the FITC channel.



Detection of change in Mitochondrial Transmembrane Potential by Mitochondrial Capture Reagent I/MitoCapture and fluorescence Microscopy: The photograph shows examples of cells scored for intact and apoptotic cells using Mitochondrial Capture Reagent I/MitoCapture probe. The apoptotic cell (left panel) shows diffused green fluorescence, whereas the normal control cell (right panel) shows punctuate red fluorescence.

6. Troubleshooting

Problem	Reason	Solution
Lower signal levels	Washing cells with PBS before/after fixation (adherent cells)	Always use binding buffer for washing cells
	Cells did not initiate apoptosis	Determine the time-point for initiation of apoptosis after induction (time-course experiment)
	Very few cells used for analysis	Refer to data sheet for appropriate cell number
	Incorrect setting of the equipment used to read samples	Refer to datasheet and use the recommended filter setting

	Use of expired kit or improperly stored reagents	Always check the expiry date and store the components appropriately
Problem	Reason	Solution
Erratic results	Uneven number of cells seeded	Seed only healthy cells (correct passage number)
	Adherent cells dislodged at the time of experiment	Perform experiment gently and in duplicates or triplicates for each treatment
	Incorrect incubation times or temperatures	Refer to datasheet & verify correct incubation times and temperatures
	Incorrect volumes used	Use calibrated pipettes and aliquot correctly

Problem	Reason	Solution
High Background	Cell density is higher than recommended	Refer to datasheet and use the suggested cell number
	Increased volumes of components added	Use calibrated pipettes accurately
	Incubation of cell samples for extended periods	Refer to datasheets and incubate for exact times
	Use of extremely confluent cells	Perform assay when cells are at 80-95% confluency
	Contaminated cells	Check for bacteria/ yeast/ mycoplasma contamination

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “contact us” on www.abcam.com for the phone number for your region).

UK, EU and ROW

Email: technical@abcam.com

Tel: +44 (0)1223 696000

www.abcam.com

US, Canada and Latin America

Email: us.technical@abcam.com

Tel: 888-77-ABCAM (22226)

www.abcam.com

China and Asia Pacific

Email: hk.technical@abcam.com

Tel: 108008523689 (中國聯通)

www.abcam.cn

Japan

Email: technical@abcam.co.jp

Tel: +81-(0)3-6231-0940

www.abcam.co.jp