

## Cell Meter™ Mitochondrial Membrane Potential Assay Kit

### \*Orange Fluorescence Optimized for Flow Cytometry\*

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 22804 (100 assays)	Keep in freezer and avoid exposure to light	Flow cytometer

### Introduction

Our Cell Meter™ assay kits are a set of tools for monitoring cell viability. There are a variety of parameters that can be used. This particular kit is designed to detect cell apoptosis by measuring the loss of the mitochondrial membrane potential (MMP). The collapse of mitochondrial membrane potential coincides with the opening of the mitochondrial permeability transition pores, leading to the release of cytochrome C into the cytosol, which in turn triggers other downstream events in the apoptotic cascade.

Our Cell Meter™ Orange Mitochondrial Membrane Potential Assay Kit provides all the essential components with an optimized assay method. This fluorimetric assay uses our proprietary cationic MitoTell™ Orange for the detection of apoptosis in cells with the loss of mitochondrial membrane potential. In normal cells, the red fluorescence intensity is increased when MitoTell™ Orange is accumulated in the mitochondria. However, in apoptotic cells, the fluorescence intensity of MitoTell™ Orange is decreased following the collapse of MMP. Cells stained with MitoTell™ Orange can be visualized with a flow cytometer at 488 nm excitation with red emission (FL2 channel). The kit can be used together with other reagents, such as Cell Meter™ Phosphatidylserine Apoptosis Assay Kit (22835) for multi-parametric study of cell vitality and apoptosis. The kit is optimized for screening apoptosis activators and inhibitors with a flow cytometer.

### Kit Components

Components	Amount
Component A: 500X MitoTell™ Orange in DMSO	1 vial (200 µL)
Component B: Assay Buffer	1 bottle (100 mL)

### Assay Protocol for Flow Cytometer

#### Brief Summary

**Prepare cells with test compounds at the density of  $5 \times 10^5$  to  $1 \times 10^6$  cells/mL → Add 2 µL of 500X MitoTell™ Orange into 1 mL of cell solution → Incubate at room temperature for 15-30 minutes  
→ Pellet the cells, and resuspend the cells in 1 mL of growth medium  
→ Analyze cells by using a flow cytometer (FL2 channel)**

*Note: Thaw all the kit components at room temperature before use.*

- For each sample, prepare cells in 1 mL of warm medium or buffer of your choice at the density of  $5 \times 10^5$  to  $1 \times 10^6$  cells/mL.  
*Note: Each cell line should be evaluated on an individual basis to determine the optimal cell density for apoptosis induction.*
- Treat cells with test compounds for a desired period of time to induce apoptosis, and set up parallel control experiments.  
For Negative Control: Treat cells with vehicle only.  
For Positive Control: Treat cells with FCCP or CCCP at 5-50 µM in a 37 °C, 5% CO<sub>2</sub> incubator for 15 to 30 minutes.  
*Note: CCCP or FCCP can be added simultaneously with MitoTell™ Orange (See Step 3). To get the best result, titration of the CCCP or FCCP may be required for each individual cell line.*

3. Add 2  $\mu$ L of 500X MitoTell™ Orange (Component A) into the treated cells (from Step 2), and incubate the cells in a 37 °C, 5% CO<sub>2</sub> incubator for 15 to 30 minutes.

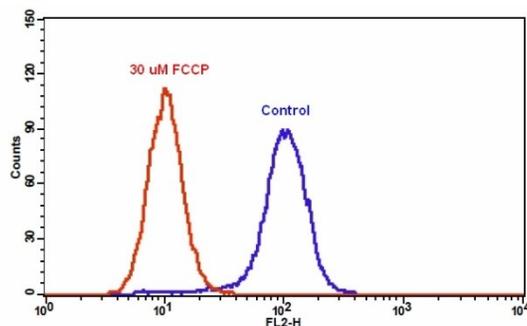
*Note: For adherent cells, gently lift the cells with 0.5 mM EDTA to keep the cells intact and wash the cells once with serum-containing media prior to the incubation with MitoTell™ Orange dye-loading solution.*

4. Centrifuge the cells at 1000 rpm for 4 minutes, and then re-suspend cells in 1 mL of Assay Buffer (Component B) or buffer of your choice.

5. Monitor the fluorescence intensity by using a flow cytometer in the FL 2 channel (Ex/Em = 540/590 nm). Gate on the cells of interest, excluding debris.

### Data Analysis

In live non-apoptotic cells, the orange fluorescence intensity is increased when MitoTell™ Orange is accumulated in the mitochondria. In apoptotic and dead cells, MitoTell™ Orange stain intensity is decreased following the collapse of MMP.



**Figure 7.43.** The decrease in fluorescence intensity of MitoTell™ Orange with the addition of FCCP in Jurkat cells. Jurkat cells were loaded with MitoTell™ Orange alone (Blue) or in the presence of 30  $\mu$ M FCCP (Red) for 15 minutes. The fluorescence intensity of MitoTell™ Orange was measured with a FACSCalibur (Becton Dickinson, San Jose, CA) flow cytometer using FL2 channel.

**Warning:** This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.