

# Mitochondrial Membrane Potential by Object Spot Counting

Using Gen5 to analyze mitochondrial  
membrane potential



## Authors

Sarah Beckman, PhD  
Agilent Technologies, Inc.  
Winooski, VT, USA

Nadia A. Rana, PhD  
Enzo Life Sciences, Inc.  
Farmingdale, NY, USA

## Abstract

Mitochondria are critical organelles involved in the maintenance of viability and vitality. However, mitochondria are also often involved in cell death and disease. These diverse functions of mitochondria are all at some point dependent on the mitochondrial membrane potential (MMP). Analysis of membrane potential is extremely valuable for obtaining insights into both basic energy metabolism and its dysfunction. This application note describes the use of a cationic MMP-sensitive fluorescent probe in combination with object-based spot counting analysis to assess the effects of carbonyl cyanide 3-chlorophenylhydrazone (CCCP) and ethanol (EtOH) on MMP. It was found that CCCP depolarized MMP in a concentration-dependent manner. In contrast, EtOH first hyperpolarized the MMP, and subsequently depolarized at concentrations above 3%.

## Introduction

The mitochondrion is a highly dynamic organelle, often called the "powerhouse of the cell" for its ability to produce cellular energy in an efficient manner. Mitochondria are the primary manufacturers of ATP, but they also regulate iron homeostasis and the production of free radicals.<sup>1</sup> Mitochondria have a duality of function in that they are involved in the maintenance of viability and vitality, but also play a role in the regulation of apoptotic cell death. Studies have demonstrated that metabolic control through mitochondria is not only related to cell fate, but also plays an important role in differentiation.<sup>2</sup> These diverse functions of mitochondria are all at some point dependent on the MMP.

Mitochondrial respiration generates an electrochemical gradient of protons made up mostly of a negative electrical potential difference across the mitochondrial inner membrane.<sup>3</sup> During mitochondrial oxidative phosphorylation, the transfer of electrons through electron transport chain (ETC) complexes I through IV in the inner mitochondrial membrane provides the energy to drive protons against their concentration gradient across the inner mitochondrial membrane (out of the mitochondrial cytoplasm). The result of this process is an accumulation of H<sup>+</sup> outside the membrane, which then flows back into the mitochondria through Complex V, producing ATP.<sup>4</sup> This accumulation of H<sup>+</sup> results in an electrochemical gradient, otherwise known as MMP.

Mitochondria are present in most of the cells in a living organism, and as such, they are implicated in a wide variety of diseases. Defects in the transfer of electrons across the mitochondrial membrane can cause electrons to accumulate on the ETC complexes and enhance reactive oxygen species (ROS) production. This accumulation increases the potential for electrons to bind with free oxygen species and contributes to many pathological conditions including degenerative diseases, cancer, and aging.<sup>5</sup> Disruption of MMP is one of the earliest intracellular events to occur following induction of apoptosis. In mammalian cells, three responses of mitochondria following a death signal have been noted: a transient hyperpolarization of MMP, a subsequent substantial depolarization of MMP, and, in selected settings, the release of cytochrome c.<sup>6,7</sup>

Cell-based assays for the analysis of MMP are extremely valuable to obtain insights into both cell disease and viability. Evaluating the functional status of mitochondria is critical to elucidating the role of mitochondrial activity in drug-induced toxicity, apoptosis, stem cells, and other cellular and biochemical processes. Quantitative microscopy of the intracellular distribution of membrane-permeant

cationic fluorophores provides a means to measure MMP in live-cultured cells.<sup>3</sup> This application note describes the use of a fluorescent probe in combination with object-based spot counting analysis to assess the effects of carbonyl cyanide 3-chlorophenylhydrazone (CCCP) and ethanol on MMP. The object spot counting capability of Agilent BioTek Gen5 3.03 software was used to perform an analysis to accurately and efficiently determine the number of MITO-ID MP-positive aggregates per cell.

## Materials and methods

### Cell culture

HeLa cells were grown in Advanced Dulbecco's Modified Eagle's Medium (DMEM) (Gibco, Grand Island, NY) with 10% FBS (Gibco) and 1x PennStrep-Glutamine (Cellgro, Manassas, VA). Cells were seeded into black-sided clear-bottom 96-well microplates (Corning, Corning, NY) at 20,000 cells per well.

### MITO-ID Membrane Potential Detection Kit

A MITO-ID Membrane Potential Detection Kit (donated by Enzo Life Sciences, Farmingdale, NY) was used to assess the change in MMP in perturbed HeLa cells. The probe is a cationic carbocyanine dye with a polychromatic fluorescent emission in which the dye fluoresces green or red, depending upon MMP. In normal healthy cells with an established proton gradient across the mitochondrial membrane, the cationic probe is largely driven into the organelle where it exists as red fluorescent aggregates; yet a portion of the dye remains as a green fluorescent monomer in the cytosol. The relative proportion of these localizations, aggregations, and resultant fluorescence is dictated by the magnitude of the MMP. Hyperpolarized cells tend to exhibit more red fluorescence. As the mitochondrial membrane depolarizes (MMP reduces), the cationic probe's tendency to follow the electrical gradient is reduced and preferentially exists as a green-fluorescent monomer in the cytosol.

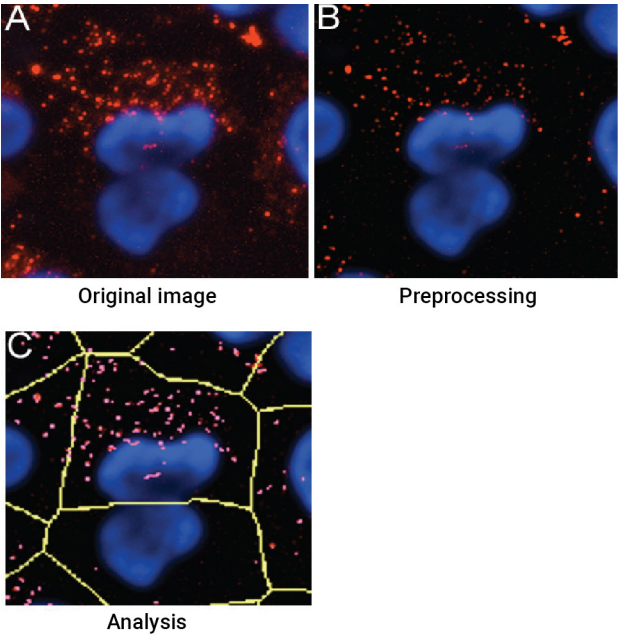
Before MITO-ID Membrane Potential (MP) detection, HeLa cells were treated either with 0.5 to 16  $\mu$ M CCCP for 30 minutes or 0.375% to 6% ethanol (EtOH) for six hours. Following treatment, cells were washed 2x with 100  $\mu$ L of 1x assay solution prepared according to the manufacturer's protocol. Next, the assay solution was replaced with 100  $\mu$ L of detection reagent (1 mL of assay solution + 2  $\mu$ L of Hoechst + 15  $\mu$ L of MITO-ID MP Detection Reagent) for 15 minutes at room temperature in the dark. The detection reagent was prepared immediately before use. Finally, the cells were washed once with assay solution and were imaged immediately.

Cell imaging

Images were acquired using a 20x objective on an Agilent BioTek Lionheart FX automated microscope configured with DAPI, GFP, and Texas Red light cubes. The DAPI light cube was configured with a 377/50 excitation filter and a 447/60 emission filter. The GFP light cube used a 469/35 excitation filter and a 525/39 emission filter. The Texas Red light cube used a 585/29 excitation filter and a 624/40 emission filter. Images were taken by automatically focusing on the DAPI channel, with an offset of 10 to 12 on the GFP and Texas Red channels.

Image analysis

Object preprocessing was used to ensure the best possible detection of nuclei and the best separation between individual aggregates. Imaging preprocessing parameters are described in detail in Table 1. The Texas Red channel of all the images was preprocessed with a 1 µm rolling ball to obtain the best separation between individual aggregates (Figure 1).



**Figure 1.** MITO-ID MP spot counting workflow. (A) Original image. (B) Preprocessed image. (C) Object masks highlighting cell area in yellow and spots in pink.

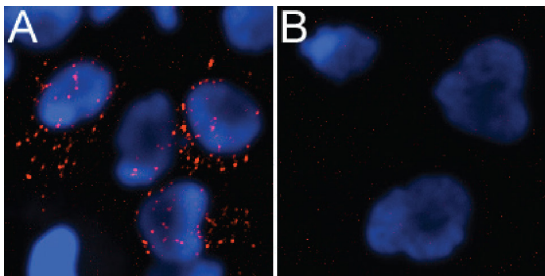
Automatic preprocessing parameters were applied to the DAPI channel to facilitate masking of the nuclei. Cell object counting analysis was then performed on the transformed DAPI channel to highlight each individual cell. Object spot counting was performed on the Texas Red channel to determine the number of MITO-ID MP aggregates per nuclei according to the parameters outlined below in Table 1.

**Table 1.** Agilent BioTek Gen5 microplate reader and imager software settings. Image analysis parameters for generating a cellular mask in the DAPI channel and an object mask in the Texas Red channel to count MITO-ID MP-positive aggregates.

| Imaging Preprocessing                        |                        |
|--|------------------------|
| Image Set                                    | DAPI                   |
| Background                                   | Dark                   |
| Rolling Bar Diameter                         | Auto                   |
| Image Smoothing Strength                     | 0                      |
| Image Set                                    | Texas Red              |
| Background                                   | Dark                   |
| Rolling Bar Diameter                         | 1 µm                   |
| Priority                                     | Fast speed             |
| Image Smoothing Strength                     | 0                      |
| Cellular Analysis                            |                        |
| Detection Channel:<br>Primary Mask and Count | Tsf[DAPI 377,447]      |
| Threshold                                    | 5,000                  |
| Secondary Mask                               | Tsf[Texas Red 586,647] |
| Measure within a Secondary Mask              | Checked                |
| Expand Primary Mask                          | 30 µm                  |
| Threshold                                    | Unchecked              |
| Count Spots                                  | Checked                |
| Size   | 0.5 to 5 µm            |
| Advanced Options                             | Count spots options    |
| Rolling Ball Size                            | Smaller                |
| Threshold                                    | 1,000                  |

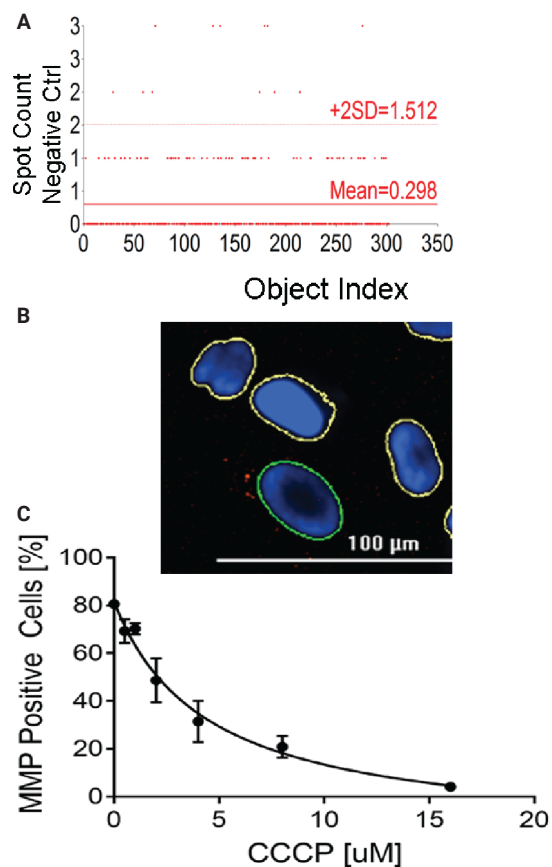
## Results and discussion

HeLa cells were treated with 0.5 to 16  $\mu\text{M}$  CCCP for 30 minutes to determine the effect of increasing concentrations of CCCP on the number of MITO-ID MP aggregates per cell. CCCP is a proton ionophore and uncoupler of oxidative phosphorylation in mitochondria. CCCP leads to a dissipation of the electrochemical potential across the inner mitochondrial membrane caused by its ability to translocate protons across membranes.<sup>8</sup> It was used in this study for depolarizing mitochondrial membranes. In this assay, decreasing red MITO-ID MP fluorescence aggregates indicate decreasing MMP as the red MITO-ID MP aggregates leave the mitochondria during depolarization. There is a decrease in the number of MMP-positive aggregates as a result of increasing CCCP concentration (Figure 2).



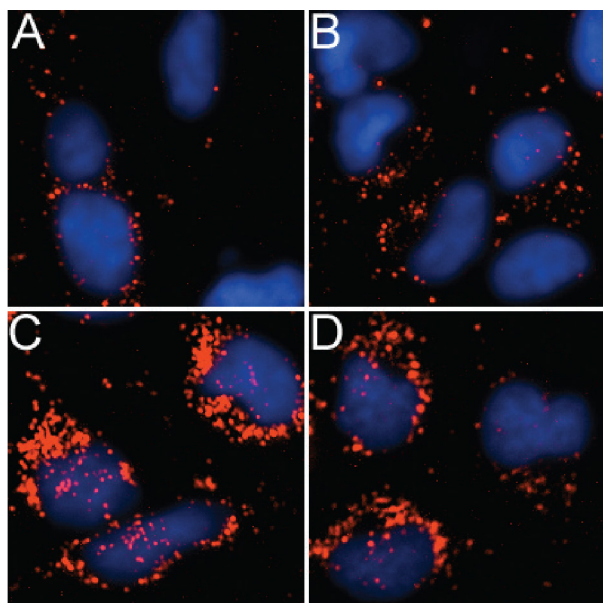
**Figure 2.** MMP-positive aggregates decrease after treatment with CCCP. (A) 0  $\mu\text{M}$  CCCP. (B) 16  $\mu\text{M}$  CCCP.

To determine the percent responders, or percent MMP-positive cells, the scatter plot function in Gen5 3.03 was used. Anything above the mean plus two standard deviations (SD) is considered a "responder". The spot count from the negative control sample was used to define what constitutes a responder. Anything above 1.5 spots per cell was determined to be a responder (Figure 3A). Figure 3B shows a responder cell highlighted in green and nonresponders highlighted in yellow. The percent of MMP-positive cells was plotted against CCCP concentration and, as expected, a decrease in percent MMP-positive cells was observed with an increase in CCCP concentration (Figure 3C).



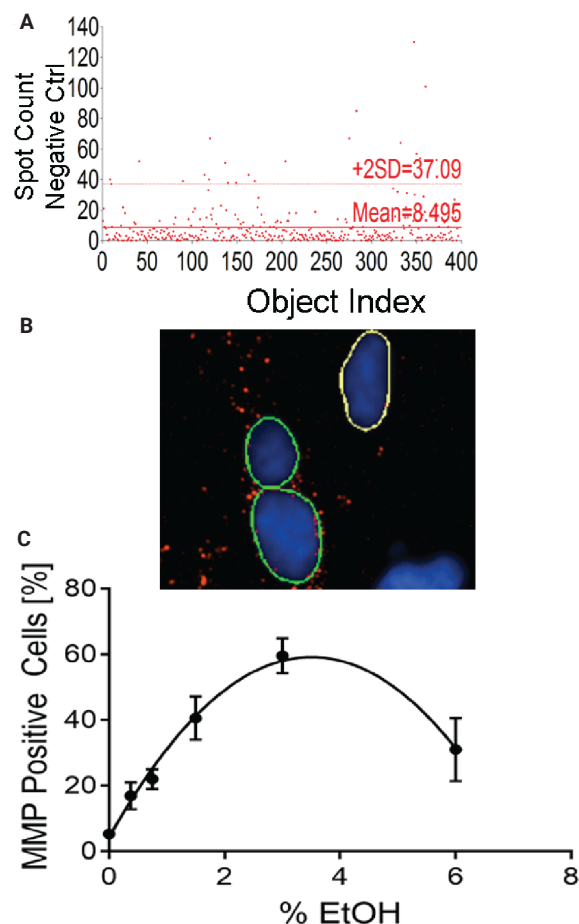
**Figure 3.** Percent MMP-positive cells following treatment with CCCP. (A) Scatter plot showing mean + 2 SD = 1.5 spots per cell. (B) Nonresponders shown in yellow, responders shown in green. (C) Percent responders decrease in a concentration-dependent manner.

HeLa cells were treated with 0.375 to 6% EtOH for 6 hours to determine the effect of increasing concentrations of EtOH on the number of MMP-positive aggregates per cell. Changes in MITO-ID MP fluorescence indicated an initial hyperpolarization of the mitochondrial membrane followed by a decrease in polarization with the highest concentration of EtOH (Figure 4). This is consistent with previous studies where ethanol at lower concentrations caused hyperpolarization of the mitochondrial membrane in ARPE-19 cells, and higher concentrations of ethanol resulted in membrane depolarization.<sup>9</sup>



**Figure 4.** MMP-positive aggregates in HeLa cells after treatment with EtOH. (A) 0  $\mu$ M EtOH, (B) 0.75% EtOH, (C) 3% EtOH, (D) 6% EtOH.

The spot count from the negative control sample of this experiment was again used to decide what constitutes a responder. It was determined that anything above 37 spots per cell was a responder (Figure 5A). Figure 5B shows responder cells highlighted in green and a nonresponder highlighted in yellow. The percent of MMP-positive cells was plotted against EtOH treatment and see an increase in percent MMP-positive cells up to 3% EtOH, then a decrease in MMP-positive cells at treatments above 3% (Figure 5C).



**Figure 5.** Percent MMP-positive cells following treatment with EtOH. (A) Scatter plot showing mean + 2 SD = 37 spots per cell. (B) Nonresponders are shown in yellow, responders are shown in green. (C) Percent responders show an initial increase with EtOH treatment followed by a decrease at concentrations above 3%.

## Conclusion

This study used MITO-ID Membrane Potential Dye in combination with an Agilent BioTek Lionheart FX automated microscope and Agilent BioTek Gen5 microplate reader and imager software to assess the effects of CCCP and EtOH treatment on the MMP of HeLa cells. MMP was quantified by the analysis of MITO-ID MP-positive aggregates in cells using the Gen5 3.03 object spot counting feature. Spot counting allows reliable and accurate measurements of object-level MITO-ID MP aggregate analysis in live-cultured cells.

## References

1. Rousselle, C. *et al.* Cytometric Assessment of Mitochondria Using Fluorescent Probes. *Cytometry Part A* **2011**, 79, 405–425.
2. Pérez, M. J. J.; Quintanilla, R. A. Development or Disease: Duality of the Mitochondrial Permeability Transition Pore. *Dev. Biol.* **2017**, 426, 1–7.
3. Lemasters, J. J.; Ramshesh, V. K. Imaging of Mitochondrial Polarization and Depolarization with Cationic Fluorophores. *Methods Cell Biol.* **2007**, 80, 283–95.
4. Perry, S. W. *et al.* Mitochondrial Membrane Potential Probes and the Proton Gradient: a Practical Usage Guide. *BioTechniques* **2011**, 50, 98–115.
5. Johannsen, D. L.; Ravussin, E. The Role of Mitochondria in Health and Disease. *Curr. Opin. Pharmacol.* **2009**, 9, 780–6.
6. Gross, A. *et al.* Biochemical and Genetic Analysis of the Mitochondrial Response of Yeast to BAX and BCL-X(L). *Mol. Cell. Biol.* **2000**, 20, 3125–36.
7. Farrace, G. M.; Piredda, L.; Matarrese, P. Transglutaminase Overexpression Sensitizes Neuronal Cell Lines to Apoptosis by Increasing Mitochondrial Membrane Potential and Cellular Oxidative Stress. *Journal of Neurochemistry* **2002**. doi:10.1046/j.1471-4159.2002.00898.x
8. Martin, J.; Mahlke, K.; Pfanner, N. Role of an Energized Inner Membrane in Mitochondrial Protein Import. Delta Psi Drives the Movement of Presequences. *Journal of Biological Chemistry* **1991**. at <<http://www.jbc.org/content/266/27/18051.short>>
9. Bonet-Ponce, L.; Saez-Atienzar, S.; da Casa, C. On the Mechanism Underlying Ethanol-Induced Mitochondrial Dynamic Disruption and Autophagy Response. *Biochimica et Biophysica Acta* **2015**.

[www.agilent.com/lifesciences/biotek](http://www.agilent.com/lifesciences/biotek)

For Research Use Only. Not for use in diagnostic procedures.

RA44412.4026157407

This information is subject to change without notice.

© Agilent Technologies, Inc. 2017, 2022  
Printed in the USA, May 12, 2022  
5994-3386EN