

MT-1 MitoMP Detection Kit

Technical Manual

Technical Manual (Japanese version) is available at <http://www.dojindo.co.jp/manual/MT13.pdf>

General Information

Mitochondria is an important organelle that uses oxygen to synthesize ATP, producing the necessary energy for living cells to thrive¹⁾. Decreased mitochondrial activity and mitochondrial dysfunction are associated with cancer, aging, and neurodegenerative diseases such as Alzheimer's and Parkinson's disease^{2,3)}. Therefore, mitochondrial membrane potential (MMP) has been widely studied as a promising target for mitochondria-related diseases.

JC-1 dye, tetramethylrhodamine ethyl ester (TMRE), and tetramethylrhodamine methyl ester (TMRM) are widely used to monitor MMP, however, these dyes have some limitations, such as low photostability and poor retention after aldehyde fixation. These limitations result in poor reproducibility of experiments.

Dojindo's MT-1 MitoMP Detection Kit overcomes these limitations. This kit can monitor MMP with excellent reproducibility even after aldehyde fixation. Furthermore, the MT-1 Dye in this kit is extremely photostable and more sensitive than JC-1. In addition, the Imaging Buffer included in this kit minimizes background fluorescence and maintains cell vitality while the assay is being performed.

Contents

MT-1 Dye 20 μ l x 3
Imaging Buffer (10x) 11 ml x 1

Storage Condition

Store at -20°C, protected from light and moisture.

Required Equipment and Materials

- Growth medium or HBSS
- Micropipettes (100–1000 μ l, 0.5–10 μ l)
- Microtubes

Preparation of Solutions

Preparation of Imaging Buffer solution

Dilute the 10x Imaging Buffer (1:10) in double-deionized water.

***Use the Imaging Buffer solution within the same day.**

Preparation of MT-1 working solution

Dilute the MT-1 Dye (1:1000) in the cell culture medium.

***Use the MT-1 working solution within the same day.**

Table 1 Required amount of Imaging Buffer solution by vessel type

Vessel	35-mm dish	ibidi 8-well plate
Appropriate amount	2 ml/dish	200 μ l/well
Vessel	96-well plate	Sample tube (Flow Cytometry)
Appropriate amount	100 μ l/well	0.5 ml/sample

General protocol

1. Inoculate cells into a dish or chamber slide and incubate in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
2. Add an appropriate volume of MT-1 working solution to cells.
3. Incubate the cells for 30 minutes in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
4. Discard the supernatant and wash the cells with medium or HBSS twice.
5. Induce a change in MMP.
6. Discard the supernatant and wash the cells with medium or HBSS twice.
7. Add Imaging Buffer solution and observe the cells under a fluorescence microscope.

Usage Examples

Fluorescence microscopic detection of MMP in HeLa cells treated with carbonyl cyanide-p-trifluoromethoxyphenyl hydrazone (FCCP)

1. HeLa cells (2.4x10⁵ cells/ml, 100 μ l) in MEM (supplemented with 10% fetal bovine serum, and 1% penicillin-streptomycin) were seeded in a 96-well clear-bottom black plate and cultured overnight in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
2. After the supernatant (100 μ l) was removed, the MT-1 working solution (1:1000 MT-1 diluted in MEM, 100 μ l) was added. The cells were cultured for 30 minutes in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
3. The cells were washed with 100 μ l of HBSS twice.
4. FCCP in MEM (0, 1, or 5 μ mol/l, 100 μ l) was added to each well, and then the cells were cultured for 30 minutes in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
5. The cells were washed with 100 μ l of HBSS twice.
6. Imaging Buffer solution (100 μ l) was added, and the cells were observed under a fluorescence microscope.

***The imaging data were quantified using Macro Cell Count (KEYENCE).**

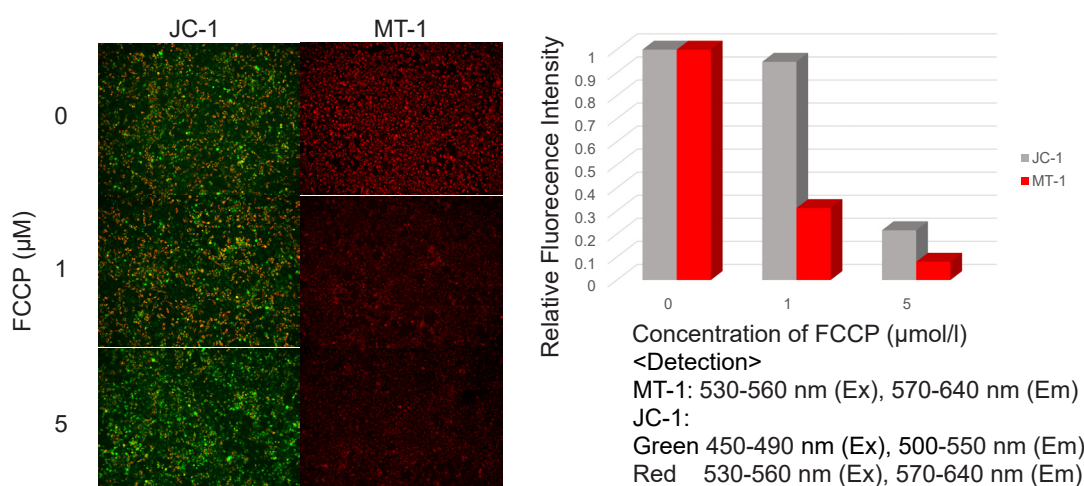


Figure 1. Fluorescence imaging of mitochondrial membrane potential in HeLa cells.

1. HL60 cells (1.0×10^6 cells/ml, 1 ml) in RPMI (10% fetal bovine serum, 1% penicillin-streptomycin) were transferred to a 1.5 ml tube.
2. The cells suspension was centrifuged at 200xg for 3 minutes, and then the medium was removed and the MT-1 working solution (1:1000 MT-1 diluted in RPMI, 1 ml) was added. Next, the cells were incubated for 30 minutes in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
3. The cells suspension was centrifuged at 200xg for 3 minutes, and then the cells were washed with 100 µl of HBSS twice.
4. Etoposide in RPMI (0 or 50 µmol/l, 1 ml) was added, and cells were incubated for 24 hours in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
5. The sample (100 µl) was transferred to another 1.5 ml tube and 5 µl of FITC Annexin V (Becton Dickinson, Cat. No. 51-65874X) was added, followed by incubation for 15 minutes in an incubator set at 37 °C and equilibrated with 95% air and 5% CO₂.
6. Imaging Buffer solution (400 µl) was added, and the cells were analyzed using a flow cytometer.

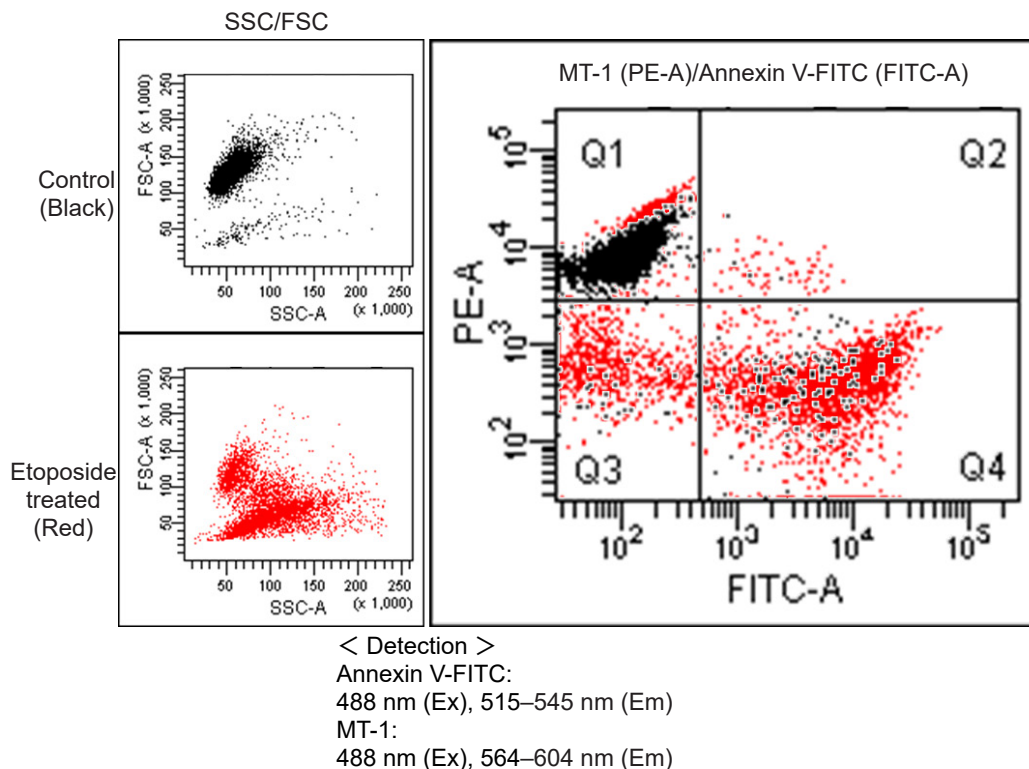


Figure 2. Flow cytometric analysis of MMP in HL60 cells.

References

- 1) Ferri, K. F. et al., *J. Exp. Med.*, **2000**, 192, 1081–1092.
- 2) Matsuda, N. et al., *J. Cell Biol.*, **2010**, 189, 211.
- 3) Wang, J. L. et al., *PNAS*, **2000**, 97, 7124–7129.

If you need more information, please contact Dojindo technical service.

Dojindo Molecular Technologies, Inc.

30 West Gude Dr., Suite 260, Rockville, MD 20850, USA

Toll free: 1-877-987-2667 Phone: 301-987-2667 Fax: 301-987-2687

E-mail: info@dojindo.com Web: www.dojindo.com

Dojindo Molecular Technologies, Inc.

Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>

Dojindo EU GmbH

Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>

Dojindo China Co., Ltd

Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>