

TMRE-Mitochondrial Membrane Potential Assay Kit ab113852

★★★★★ [4 Abreviews](#) [166 References](#) [画像数 6](#)

医薬用外劇物

製品の概要

製品名	TMRE-Mitochondrial Membrane Potential Assay Kit
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Cell-based (qualitative)
全工程の試験時間	0h 35m
製品の概要	Store kit at 4°C in the dark immediately upon receipt.

TMRE-Mitochondrial Membrane Potential Assay Kit ab113852 is used for quantifying changes in mitochondrial membrane potential in live cells by flow cytometry, microplate spectrophotometry and fluorescent microscopy.

TMRE (tetramethylrhodamine, ethyl ester) is used to label active mitochondria. TMRE is a cell permeant, positively-charged, red-orange dye that readily accumulates in active mitochondria due to their relative negative charge. Depolarized or inactive mitochondria have decreased membrane potential and fail to sequester TMRE. NB: TMRE is also available as free molecule as [ab274305](#) (Tetramethylrhodamine ethyl ester).

The TMRE protocol also uses FCCP (carbonyl cyanide 4-(trifluoromethoxy) phenylhydrazone), which is a ionophore uncoupler of oxidative phosphorylation. Treating cells with FCCP eliminates mitochondrial membrane potential and TMRE staining. TMRE is suitable for the labeling of mitochondria in live cells and is not compatible with fixation.

TMRE protocol summary:

- add FCCP to appropriate control cell samples and incubate for 10 min
- incubate with TMRE for 15-30 min, pellet (suspension cells) / remove media (adherent cells) and wash with PBS / 0.2% BSA
- analyze with micro-plate reader at Ex/Em 549/575 nm, flow cytometer using 488nm laser for excitation and at emission 575 nm, or fluorescent microscope.

特記事項

TMRE is only suitable for use with live (not fixed) cells.

Related assays

Review the [cell health assay guide](#) to learn about kits to perform a [cell viability assay](#), [cytotoxicity assay](#) and [cell proliferation assay](#).

Review the [metabolism assay guide](#) to learn about assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also about how to assay metabolic function in live cells using your plate reader.

How other researchers have used TMRE assay kit ab113852

Maiti AK et al. of the University of Gothenburg, Sweden, used TMRE assay kit ab113852 to identify that *C. rodentium* infection in mice* reduced mitochondrial transmembrane potential. In a subsequent paper, they identified that this effect was reduced by treatment with vasoactive intestinal peptide (VIP). *a model for enteropathogenic *E. coli*

試験プラットフォーム

Microplate reader, Fluor. microscope, Flow cyt.

製品の特性

保存方法

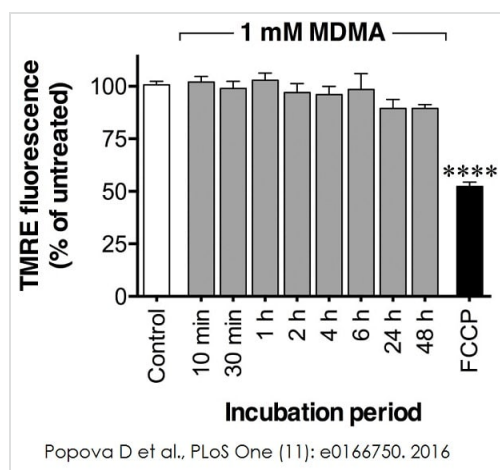
Store at +4°C. Please refer to protocols.

内容	200 tests
1mM TMRE (in DMSO)	1 x 40µl
50mM FCCP (in DMSO)	1 x 10µl

関連性

Mitochondrial Membrane Potential is an important parameter of mitochondrial function used as an indicator of cell death. The collapse of the 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.

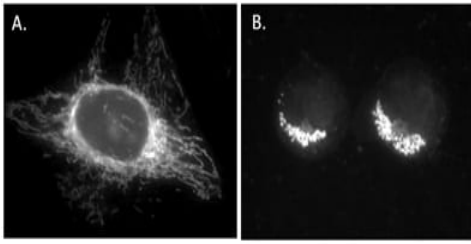
画像



P19 neurons (750 cells/mm²) were exposed to MDMA on days 7–9 in serum-free medium for 10 min up to 48 hours. The positive control FCCP (carbonyl cyanide-*p*-trifluoromethoxyphenylhydrazone), an uncoupler of mitochondrial oxidative phosphorylation, was applied at the concentration of 5 µM for 10 min. The cells were incubated with 500 µM TMRE for 30–45 min at 37°C, 5% CO₂, followed by washing once with 100 µl of HBSS containing 0.2% bovine serum albumin. A volume of 200 µL of HBSS containing 0.2% bovine serum albumin was added to each well, and the fluorescence was measured with excitation/emission: 544/590 nm.

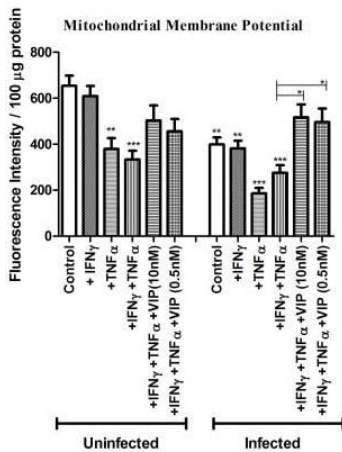
Mitochondrial membrane potential was measured using ab113852

Popova D et al., PLoS One, 11(11). Fig 5. doi: 10.1371/journal.pone.0166750 Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>



Immunocytochemistry/ Immunofluorescence - TMRE
—Mitochondrial Membrane Potential Assay Kit
(ab113852)

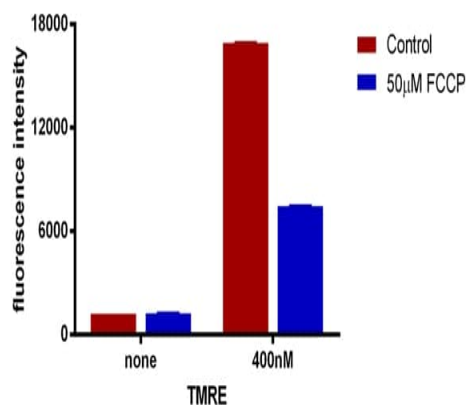
A: HeLa cells (adherent) were cultured on coverslips and stained with ab113852 (200nM TMRE) for 20 minutes in media, washed briefly with PBS and immediately imaged. **B:** Jurkat cells (suspension) were stained and washed as above and then transferred to a slide and immobilized under a coverslip for imaging.



Mitochondrial membrane potential assayed using
TMRE in the study of VIP

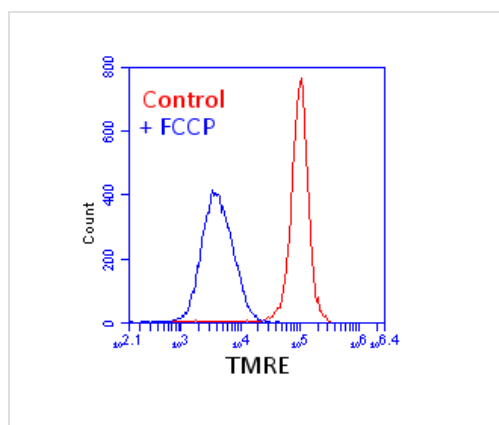
Image courtesy of Maiti AK et al. PLoS One. 2018;
13(9): e0204567. doi: 10.1371/journal.pone.0204567

Maiti AK et al (2018) used TMRE assay kit ab113852 to measure mitochondrial membrane potential in an in vitro mouse intestinal model treated with cytokines in the presence and absence of VIP (vasoactive intestinal peptide). VIP was induced by *C. rodentium* infection and cytokines.



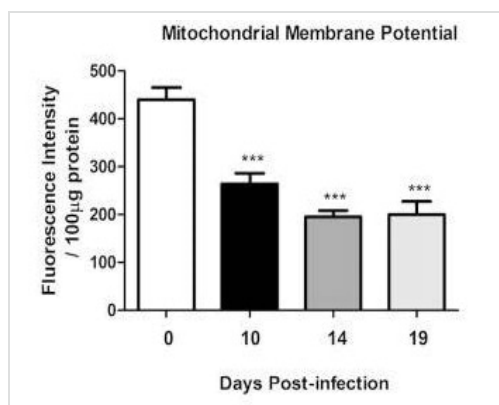
Analysis of TMRE staining using a fluorescent plate
reader and a microplate.

Chart showing mean fluorescent intensity \pm standard deviation from quadruplicate measurements of 400 nM TMRE stained Jurkat cells in a 96-well microplate \pm treatment with FCCP.



Flow cytometry histogram of Jurkat cells stained with ab113852 (100nM TMRE) with (blue) or without (red) treatment with 100µM FCCP.

Flow Cytometry - TMRE—Mitochondrial Membrane Potential Assay Kit (ab113852)



Maiti et al (2015) used TMRE assay kit ab113852 to assess mitochondrial membrane potential in murine distal colon after *C. rodentium* infection.

TMRE assay to assess mitochondrial function after *C. rodentium* infection

Image courtesy of Maiti AK et al. Sci Rep. 2015; 5: 1543. doi: 10.1038/srep15434. Reproduced under the Creative Commons License <http://creativecommons.org/licenses/by/4.0/>

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors