

ab112150

NIR Mitochondrial Membrane Potential Assay Kit (Microplate)

Instructions for Use

For measuring Mitochondrial membrane potential in cells using our proprietary fluorescence probe

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

Table of Contents

| 1. | Introduction | 3 |
|----|----------------------|----|
| 2. | Protocol Summary | 4 |
| 3. | Kit Contents | 5 |
| 4. | Storage and Handling | 5 |
| 5. | Assay Protocol | 6 |
| 6. | Data Analysis | ç |
| 7. | Troubleshooting | 10 |

1. Introduction

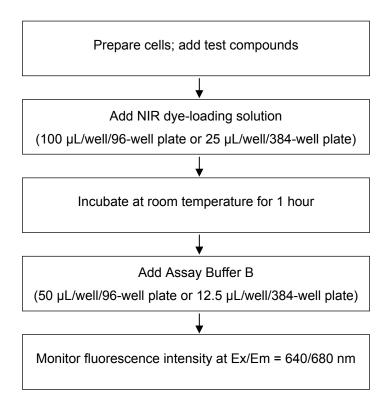
ab112150 is designed to detect cell apoptosis by measuring the loss of the mitochondrial membrane potential. The collapse of mitochondrial membrane potential coincides with the opening of the mitochondrial permeability transition pores, leading to the release of Cytochrome C into the cytosol, which in turn triggers other downstream events in the apoptotic cascade.

ab112150 NIR Mitochondria Membrane Potential Assay Kit provides all the essential components with an optimized assay method. This fluorometric assay uses our proprietary cationic a MitoNIR Dye for the detection of the mitochondrial membrane potential change in cells. In normal cells, the red fluorescence intensity is increased when MitoNIR Dye is accumulated in the mitochondria. However, in apoptotic cells, NIR stain intensity is decreased following the collapse of MMP. Cells stained with MitoNIR Dye can be monitored fluorometrically at 660-680 nm with excitation at 620-640 nm.

ab112150 can be used for screening apoptosis activators and inhibitors. The assay can be performed in a convenient 96-well and 384-well fluorescence microtiter-plate format

2. Protocol Summary

Summary for One 96-well Plate



Note: Thaw all the kit components to room temperature before starting the experiment.

3. Kit Contents

| Components | Amount |
|---------------------------------------|------------|
| Component A: 200X MitoNIR Dye in DMSO | 1 x 250 μL |
| Component B: Assay Buffer A | 1 x 50 mL |
| Component C: Assay Buffer B | 1 x 25 mL |

4. Storage and Handling

Keep at -20°C. Avoid exposure to light.

5. Assay Protocol

Note: This protocol is for one 96 - well plate.

A. Preparation of Cells

- For adherent cells: Plate cells overnight in growth medium at 20,000 to 80,000 cells/well/100 μL for a 96-well plate or 5,000 to 20,000 cells/well/25 μL for a 384-well plate.
- 2. For non-adherent cells: Centrifuge the cells from the culture medium and then suspend the cell pellets in culture medium at 100,000-200,000 cells/well/90 μL for a 96-well poly-D lysine plate or 25,000-50,000 cells/well/20 μL for a 384-well poly-D lysine plate. Centrifuge the plate at 800 rpm for 2 minutes with brake off prior to the experiments.

Note: Each cell line should be evaluated on an individual basis to determine the optimal cell density for apoptosis induction

B. Preparation of NIR Dye-loading Solution

- Thaw all the kit components at room temperature before use.
- Add 50μL of MitoNIR Dye (Component A) into 10 mL of Assay Buffer A (Component B), and mix them well.

Note: Aliquot and store the unused MitoNIR Dye (Component A) at -20°C. Avoid repeated freeze/thaw cycles.

C. Run MitoNIR Dye Assay

 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 μ M in a 37°C, 5% CO₂ incubator for 15 to 30 minutes.

Note: CCCP or FCCP can be added simultaneously with NIR Dye. To get the best result, titration of the CCCP or FCCP may be required for each individual cell line.

2. Remove the cell medium before adding NIR dye-loading solution (See Step C.3).

Note: It is important to remove the cell medium before adding NIR dye-loading solution.

3. Add 100 μ L/well/96-well plate or 25 μ L/well/384-well plate of NIR dye-loading solution (from Step B.2) into the cell plate (from Step C.2).

4. Incubate the dye-loading plate in a 37°C, 5% CO₂ incubator for 15-30 minutes, protected from light.

Note: The appropriate incubation time depends on the individual cell type and cell concentration used. Optimize the incubation time for each experiment.

 Add 50 μL/well/96-well plate or 12.5 μL/well/384-well plate of Assay Buffer B (Component C) into the dye-loaded cell plate (from Step C.4) before monitoring the fluorescence signal.

Note 1: DO NOT wash the cells after loading.

Note 2: For non-adherent cells, it is recommended to centrifuge cell plates at 800 rpm for 2 minutes with brake off after adding Assay Buffer B (Component C).

Monitor the fluorescence intensity at Ex/Em = 640/680 nm (bottom read) either using the endpoint mode or using the kinetic mode 10 to 30 minutes after Step C.5.

6. Data Analysis

In live non-apoptotic cells, the red fluorescence intensity is increased when the MitoNIR Dye is accumulated in the mitochondria. In apoptotic and dead cells, NIR stain intensity is decreased following the collapse of MMP.

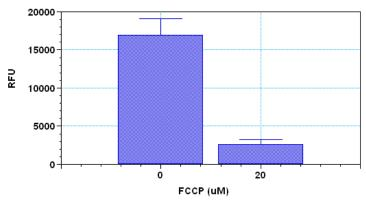


Figure 1. The decrease in NIR fluorescence with the addition of FCCP in HeLa cells. HeLa cells were dye loaded with MitoNIR Dye alone or in the presence of 20 μ M FCCP for 15 minutes. The fluorescence intensity of MitoNIR Dye was measured 30 minutes after adding Assay buffer B (Component C) with a microplate reader at Ex/Em = 640/680 nm (cut off 665 nm, bottom read).

7. Troubleshooting

| Problem | Reason | Solution |
|--------------------|------------------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Assay not working | Assay buffer at wrong temperature | Assay buffer must not be chilled - needs to be at RT |
| | Protocol step missed | Re-read and follow the protocol exactly |
| | Plate read at incorrect wavelength | Ensure you are using appropriate reader and filter settings (refer to datasheet) |
| | Unsuitable microtiter plate for assay | Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells |
| Unexpected results | Measured at wrong wavelength | Use appropriate reader and filter settings described in datasheet |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Unsuitable sample type | Use recommended samples types as listed on the datasheet |
| | Sample readings are outside linear range | Concentrate/ dilute samples to be in linear range |

| Problem | Reason | Solution |
|---------------------------------|----------------------------------------------------------------|-----------------------------------------------------------------------------------------------------------|
| Samples with | Unsuitable sample type | Refer to datasheet for details about incompatible samples |
| inconsistent readings | Samples prepared in the wrong buffer | Use the assay buffer provided (or refer to datasheet for instructions) |
| | Samples not deproteinized (if indicated on datasheet) | Use the 10kDa spin column (ab93349) or Deproteinizing sample preparation kit (ab93299) |
| | Cell/ tissue samples not sufficiently homogenized | Increase sonication time/ number of strokes with the Dounce homogenizer |
| | Too many freeze- thaw cycles | Aliquot samples to reduce the number of freeze-thaw cycles |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Samples are too old or incorrectly stored | Use freshly made samples and store at recommended temperature until use |
| Lower/ Higher readings in | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| samples and standards | Out-of-date kit or incorrectly stored reagents | Always check expiry date and store kit components as recommended on the datasheet |
| | Reagents sitting for extended periods on ice | Try to prepare a fresh reaction mix prior to each use |
| | Incorrect incubation time/ temperature | Refer to datasheet for recommended incubation time and/ or temperature |
| | Incorrect amounts used | Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume) |

| Standard curve is not linear | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
|------------------------------------|--------------------------------------------------------|--------------------------------------------------------------------------------------------------------------------------------|
| | Pipetting errors when setting up the standard curve | Try not to pipette too small volumes |
| | Incorrect pipetting when preparing the reaction mix | Always prepare a master mix |
| | Air bubbles in wells | Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates |
| | Concentration of standard stock incorrect | Recheck datasheet for recommended concentrations of standard stocks |
| | Errors in standard curve calculations | Refer to datasheet and re-check the calculations |
| | Use of other reagents than those provided with the kit | Use fresh components from the same kit |

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: 400 921 0189 / +86 21 2070 0500

www.abcam.cn

Japan

Email: technical@abcam.co.jp

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

www.abcam.co.jp

Copyright © 2012 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.