

MITO-ID[®] Green Detection Kit

Catalog #: ENZ-51022

ENZ-51022-K500 500 assays

ENZ-51022-0100 100 assays



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Introduction	2
Precautions	3
Reagents Provided and Storage	3
Additional Materials Required	4
Methods and Procedures	4
Reagent Preparation	4
Staining Live, Adherent Cells	5
Staining Live Cells Grown in Suspension	5
Staining of Aldehyde/Detergent Permeabilized Cells	6
Appendices	7
Results	7
References	8
Trouble Shooting	9
Contact Information	10

INTRODUCTION

Enzo Life Sciences' Mito-ID[®] Green Detection Kit contains a novel mitochondria-selective dye suitable for live-, detergent-permeabilized and aldehyde-fixed cell staining. Conventional fluorescent stains for mitochondria, such as JC-1 (Cat. No. ENZ-52304), rhodamine 123 (Cat. No. ENZ-52307) and tetramethylrhodamine ethyl ester (TMRE, Cat. No. ENZ-52309), are readily sequestered by actively respiring mitochondria. However, these mitochondrial stains are subsequently leached out of cells once the mitochondria's membrane potential dissipates. This characteristic severely limits their use in experiments in which cells must be treated with nonionic detergents, aldehyde fixatives or other agents that affect the energetic state of the mitochondria. Micromolar concentrations of MITO-ID[®] Green dye are sufficient for staining mammalian cells. This has been validated with a human cervical carcinoma cell line, HeLa, a human T-lymphocyte cell line, Jurkat, and human bone osteosarcoma epithelial cell line, U2OS.

One important application of MITO-ID[®] Green dye is in fluorescence co-localization imaging with red fluorescent protein (RFP)-tagged proteins. This is a powerful approach for determining the targeting of molecules to intracellular compartments, and for screening of associations and interactions between these molecules. Additionally, many organelle-targeting probes photobleach rapidly, are subject to quenching when concentrated in organelles, are highly toxic, or only transiently associate with the target organelle, requiring imaging within a minute or two of dye addition.^{3,4}

Enzo's MITO-ID[®] Green dye, a cell-permeable small organic probe molecule that spontaneously localizes to live or fixed mitochondria, was developed to overcome the above problems. The MITO-ID[®] Green dye can be readily used in combination with other common UV and visible light excitable fluorescent dyes and various fluorescent proteins in multi-color imaging and detection applications. It emits in the FITC region of the visible light spectrum, and is highly resistant to photobleaching and concentration quenching.

The MITO-ID[®] Green Detection Kit has been specifically designed for use with RFP-expressing cell lines, as well as cells expressing blue or cyan fluorescent proteins (BFPs and CFPs). Additionally, the kit is suitable for use with live or fixed cells in conjunction with fluorescent probes, such as labeled antibodies, or other fluorescent conjugates displaying similar spectral properties as

coumarin, Texas Red and cyanine 5. A nuclear counterstain, Hoechst 33342, is provided to highlight this organelle as well.

SAFETY WARNINGS & PRECAUTIONS

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.



Handle with care

- This product is for research use only and is not intended for diagnostic purposes.
- The [®] Green Detection Reagent contains DMSO which is readily absorbed through the skin. It is harmful if ingested or absorbed through the skin and may cause irritation to the eyes. Observe appropriate precautions when handling.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas. All blood components and biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.
- To avoid photobleaching, perform all manipulations in low light environments or protected from light by other means.

REAGENTS PROVIDED AND STORAGE

All reagents are shipped on dry ice. Upon receipt, the kit should be stored at $\leq -20^{\circ}\text{C}$, protected from light. When stored properly, these reagents are stable for at least twelve months. Avoid repeated freezing and thawing.

Reagents provided in the ENZ-51022-K500 kit are sufficient for approximately 500 assays and reagents provided in the ENZ-51022-0100 kit are sufficient for approximately 100 assays using either live, adherent cells or cells in suspension.



Reagents require separate storage conditions.

Reagent	ENZ-51022-K500	ENZ-51022-0100
MITO-ID [®] Green Detection Reagent	50 μl	13 μl
Hoechst 33342 Nuclear Stain	50 μl	13 μl
10X Assay Buffer	15ml	3ml

ADDITIONAL MATERIALS REQUIRED

1. Standard fluorescence microscope
2. Calibrated, adjustable precision pipetters, preferably with disposable plastic tips
3. Adjustable speed centrifuge with swinging buckets (for suspension cultures)
4. Glass microscope slides
5. Glass cover slips
6. Deionized water
7. Anhydrous DMSO (optional)
8. Growth medium (e.g., Dulbecco's Modified Eagle Medium, D-MEM)
9. Formaldehyde (optional, for fixation protocol)
10. Triton X-100 (optional, for permeabilization protocol)

METHODS AND PROCEDURES

NOTE: Allow all reagents to thaw at room temperature before starting with the procedures. Upon thawing, gently hand-mix or vortex the reagents prior to use to ensure a homogenous solution. Briefly centrifuge the vials at the time of first use, as well as for all subsequent uses, to gather the contents at the bottom of the tube.

REAGENT PREPARATION

1. 1X Assay Buffer

Allow the 10X Assay Buffer to warm to room temperature. Make sure that the reagent is free of any crystallization before dilution. Prepare enough 1X Assay Buffer for the number of samples to be assayed by diluting each milliliter (ml) of the 10X Assay Buffer with 9ml of deionized water.

2. Dual Detection Reagent

The concentration of MITO-ID[®] Green dye for optimal staining will vary depending upon the application. Suggestions are provided to use as guidelines, though some modifications may be required depending upon the particular cell type employed and other factors such as the permeability of the dye to the cells or tissues. To reduce potential artifacts from overloading of the cells, the concentration of the dye should be kept as low as possible.

Prepare sufficient amount of Dual Detection Reagent for the number of samples to be assayed as follows: For every 1ml of 1X Assay Buffer (see preparation in step 1) or cell culture medium, add 1µl of MITO-ID® Green Detection Reagent and 1µl of Hoechst 33342 Nuclear Stain.

NOTE:

- (a) *The dyes may be combined into one staining solution or each may be used separately, if desired.*
- (b) *The Hoechst 33342 Nuclear Stain can be diluted further if its staining intensity is much stronger than the green Mitochondrial stain, MITO-ID® Green.*
- (c) *When staining BFP- or CFP-expressing cells, the Hoechst 33342 Nuclear Stain should be omitted due to its spectral overlap with these fluorescent proteins.*

STAINING LIVE, ADHERENT CELLS

1. Grow cells on cover slips inside a Petri dish filled with the appropriate culture medium. When the cells have reached the desired level of confluence, carefully remove the medium.
2. Dispense sufficient volume of Dual Detection Reagent (see Reagent Prep, section 2, page 4) to cover the monolayer cells (~100µl of labeling solution for cells grown on an 18 X 18 mm coverslip).
3. Protect samples from light and incubate for 15-30 minutes at 37°C.
4. (Optional) Wash the cells with 100µl 1X Assay Buffer. Remove excess buffer and place coverslip on slide.
5. Analyze the stained cells by wide-field fluorescence or confocal microscopy (60X magnification recommended). Use a standard FITC filter set for imaging the mitochondria. Optionally, image the nucleus using a DAPI filter set and the RFP-tagged protein using a Texas Red filter set.

STAINING LIVE CELLS GROWN IN SUSPENSION

1. Centrifuge cells for 5 minutes at 400 x g at room temperature (RT) to obtain a cell pellet.
2. Carefully remove the supernatant by aspiration and dispense sufficient volume of Dual Detection Reagent (see Reagent Prep, section 2, page 4) to cover the dispersed cell pellet.
3. Protect samples from light and incubate for 15 to 30 minutes at 37°C.

4. (Optional) Wash the cells with 100µl 1X Assay Buffer. Remove excess buffer. Resuspend cells in 30µl 1X Assay Buffer, then apply the cells to a glass slide and overlay with a coverslip.
5. Analyze the stained cells by wide-field fluorescence or confocal microscopy (60X magnification recommended). Use a standard FITC filter set for imaging the mitochondria. Optionally, image the nucleus using a DAPI filter set and the RFP-tagged protein using a Texas Red filter set.

STAINING OF ALDEHYDE-FIXED AND DETERGENT PERMEABILIZED CELLS

The MITO-ID® Green dye is capable of staining already fixed and permeabilized cells.

1. Growth of cells should be performed as described in previous two sections.
2. Carefully remove the growth medium, or 1X Assay Buffer covering the cells, and replace it with freshly prepared medium or buffer containing 3.7% formaldehyde.
3. Incubate the cells at 37°C for 15 minutes.
4. After fixation, wash the cells in PBS or 1X Assay Buffer.
5. If the cells are to be subsequently labeled with an antibody, a permeabilization step is usually required to enhance the antigen's accessibility. Incubate the fixed cells in PBS or 1X Assay Buffer containing 0.1% Triton X-100 at room temperature for about one minute.
6. Following permeabilization, rinse the cells twice in PBS or 1X Assay Buffer.
7. Perform staining as recommended for adherent or suspension cells (in previous two section) using a 500-fold dilution of the MITO-ID® Green instead of the 1000-fold dilution.

NOTE: *If desired, standard immunofluorescence staining protocols using coumarin-based, cyanine 5 or other red or blue fluorescent antibody conjugates, or equivalent, should be performed before Staining with MITO-ID® Green. Antifade compounds or mounting media may be beneficial.*

Try to view the samples as soon as possible after staining for sharper staining.

APPENDICES

FILTER SET SELECTION

The selection of optimal filter sets for a fluorescence microscopy application requires matching the optical filter specifications to the spectral characteristics of the dyes employed in the analysis. Consult the microscope or filter set manufacturer for assistance in selecting optimal filter sets for your microscope.

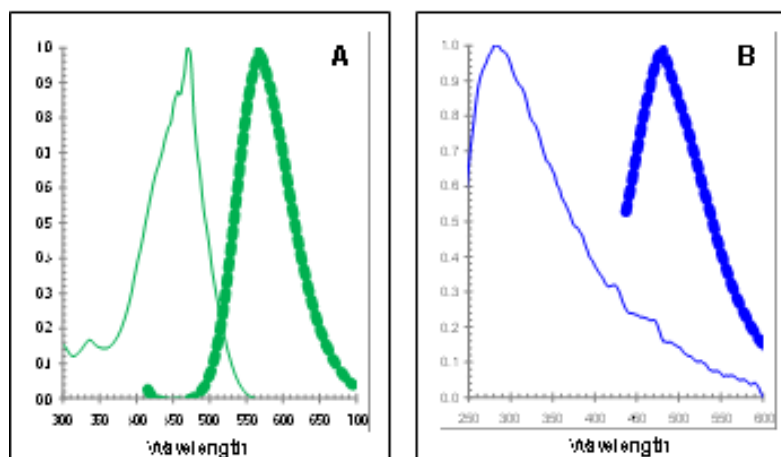


Figure 1. Absorption/excitation and fluorescence emission spectra for MITO-ID[®] Green [Ex/Em: 460/560nm] (panel A) and Hoechst 33342 [Ex/Em: 350/461nm] (panel B) dyes. All spectra were determined in 1X Assay Buffer.

RESULTS

Mitochondria are subcellular organelles found in eukaryotic cells, often representing as much as 10% of the total cell volume. Although conventional fluorescent stains for mitochondria, such as JC-1, rhodamine 123 and tetramethylrhodamine, are readily sequestered by functioning mitochondria, they are subsequently leached out of the cells once the mitochondrial membrane potential is dissipated. MITO-ID[®] Green dye accumulates in the mitochondria regardless of the mitochondrial membrane potential. The dye selectively stains mitochondria of living cells and is relatively insensitive to mitochondrial membrane potential uncouplers of phosphorylation, such as CCCP (carbonyl cyanide 3-chlorophenylhydrazone), as well as ion-channel drugs, such as valinomycin. In addition to being a live cell-permeable dye, the MITO-ID[®] Green dye is functional after cell fixation and detergent permeabilization.

MITO-ID[®] Green dye has been shown to co-localize with MITO-ID[®] Red dye. Previously, the MITO-ID[®] Red dye has been shown to co-localize with mitochondria expressing a mitochondrial-GFP tag.

REFERENCES

1. Freundt, Czapiga and Lenardo (2007) "Photoconversion of LysoTracker Red to a green fluorescent molecule" *Cell Res.* 17(11):956-958.
2. Nadrigny, Li, Kemnitz, Ropert, Koulakoff, Rudolph, Vitali, Giaume, Kirchhoff and Oheim (2007) "Systematic colocalization errors between acridine orange and EGFP in astrocyte vesicular organelles" *Biophys J.* 93(3):969-980.
3. Minamikawa, Sriratana, Williams, Bowser, Hill and Nagley (1999) Chloromethyl-X-rosamine (MitoTracker Red) photosensitises mitochondria and induces apoptosis in intact human cells. *Journal of Cell Science* 112, 2419-2430.
4. Scorrano, Petronilli, Colonna, Di Lisa and Bernardi (1999) Chloromethyltetramethylrosamine (Mitotracker Orange®) Induces the Mitochondrial Permeability Transition and Inhibits Respiratory Complex I: Implications for the mechanism of cytochrome c release" *J. Biol Chem.* 274, 35, 24657–24663.

TROUBLESHOOTING GUIDE

Problem	Potential Cause	Suggestion
Mitochondria are not sufficiently stained.	Very low concentration of MITO-ID [®] green dye was used or dye was incubated with the cells for an insufficient length of time.	Either increase the labeling concentration or increase the time allowed for the dye to accumulate in the mitochondria.
Precipitate is seen in the 10X Assay Buffer.	Precipitate forms at low temperatures.	Allow solution to warm to room temperature or 37°C, then vortex to dissolve all precipitate.
Blue nuclear counterstain is too bright compared to the green mitochondrial stain.	Different microscopes, cameras and filters may make some signals appear very bright.	Reduce the concentration of the nuclear counterstain or shorten the exposure time.
Cells do not appear healthy	Some cells require serum to remain healthy.	Add serum to stain and wash solutions. Serum does not affect staining. Normal amounts of serum added range from 2% to 10%.



Product Manual

www.enzolifesciences.com

Enabling Discovery in Life Science®

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com