



ROS-ID[®] Total ROS/Superoxide Detection Kit

**for fluorescence microscopy, flow cytometry and
microplate assay**

Instruction Manual

Catalog Number: **ENZ-51010**

200 fluorescence microscopy assays or 50 flow cytometry
assays or 2 x 96-well plates



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
Reagents Provided and Storage	3
Other Materials Needed	3
Safety Warnings and Precautions	4
Methods and Procedures	4
Reagent Preparation.....	4
Cell Preparation	6
Fluorescence/Confocal Microscopy (Adherent cells).....	6
Fluorescence/Confocal Microscopy (Suspension cells)	7
Flow Cytometry (Adherent cells)	8
Flow Cytometry (Suspension cells)	10
Fluorescence Microplate Assay (Adherent cells)	11
Fluorescence Microplate Assay (Suspension cells)	12
Appendices	14
Filter Set Selection.....	14
Setting up Optimal Exposure Time, Detection of the Dye	14
Compensation Correction for Flow Cytometry.....	15
Anticipated Results: Fluorescence Microscopy	16
Flow Cytometry Data Analysis, Anticipated Results	17
Fluorescence Microplate Assay Analysis, Results	19
References.....	21
Troubleshooting Guide	22
Contact Information.....	26

I. INTRODUCTION

Free radicals and other reactive species play influential roles in many human physiological and pathophysiological processes, including cell signaling, aging, cancer, atherosclerosis, macular degeneration, sepsis, various neurodegenerative diseases (Alzheimer's and Parkinson's disease) and diabetes. Once produced within a cell, free radicals can damage a wide variety of cellular constituents, including proteins, lipids and DNA. However, at lower concentrations these very same agents may serve as second messengers in cellular signaling. Information-rich methods are required to quantify the relative levels of various reactive species in living cells and tissues, due to the seminal role they play in physiology and pathophysiology. The ROS-ID® Total ROS/Superoxide Detection Kit enables detection of comparative levels of total ROS/RNS and also allows determination of superoxide production in live cells. Through the combination of two specific fluorescent probes, the kit provides a simple and specific assay for the real-time measurement of global levels of reactive oxygen species (ROS), peroxy nitrite and specifically superoxide in living cells.

This kit is designed to directly monitor real time reactive oxygen and/or superoxide species production in live cells using fluorescence microscopy and/or flow cytometry. The kit includes two fluorescent dye reagents as major components: Oxidative Stress Detection Reagent (Green) for ROS detection and Superoxide Detection Reagent (Orange). The non-fluorescent, cell-permeable ROS detection dye (green probe) reacts directly with a wide range of reactive species, such as hydrogen peroxide (H_2O_2), peroxy nitrite ($ONOO^-$), hydroxyl radicals (HO), nitric oxide (NO), and peroxy radical (ROO), yielding a green fluorescent product indicative of cellular production of different ROS/RNS types. However, the green probe has a low sensitivity for superoxide (O_2^-), as well as Hypochlorous acid ($HClO$) and nitric oxide (NO), relative to other free radicals. The superoxide detection dye (orange probe) is a cell-permeable probe that reacts specifically with superoxide (O_2^-), generating an orange fluorescent product. The kit is not designed to detect reactive chlorine or bromine species, as the fluorescent probes included are relatively insensitive to these analytes. Upon staining, the fluorescent products generated by the two dyes can be visualized using a

wide-field fluorescence microscope equipped with standard green (490/525 nm) and orange (550/620 nm) filter set, or cytometrically using any flow cytometer equipped with a blue (488 nm) laser.

II. REAGENTS PROVIDED AND STORAGE

All reagents are shipped on dry ice. Upon receipt, the kit should be stored at -20°C, or -80°C for long term storage. When stored properly, these reagents are stable for at least twelve months. **Avoid repeated freezing and thawing.**

Reagents provided in the kit are sufficient for at least 200 microscopy assays or 50 flow cytometry assays using live cells (adherent or in suspension)

Reagent	Quantity
Oxidative Stress Detection Reagent (Green)	300 nmol
Superoxide Detection Reagent (Orange)	300 nmol
ROS Inducer (Pyocyanin)	1 µmol
ROS Inhibitor (N-acetyl-L-cysteine)	2 x 10 mg
Wash Buffer Salts	1 pack

III. OTHER MATERIALS NEEDED

- CO₂ incubator (37°C)
- Standard fluorescence microscope or flow cytometer equipped with a blue laser (488 nm)
- Calibrated, adjustable precision pipettors, preferably with disposable plastic tips
- 5 mL round bottom polystyrene tubes for holding cells during induction of ROS/RNS (for suspension cells only) and during staining and assay procedure
- Adjustable speed centrifuge with swinging buckets
- Glass microscope slides
- Glass cover slips
- Deionized water
- Anhydrous DMF (100%)

IV. SAFETY WARNINGS AND PRECAUTIONS



Handle
with care



Protect
from light

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

V. METHODS AND PROCEDURES

NOTE: Allow all reagents to warm to room temperature before starting with the procedures. Upon thawing of solutions, 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.

A. REAGENT PREPARATIONS

Reconstitution or dilution of any and all reagents in DMSO should be avoided, as this solvent inhibits hydroxyl radical generation in cells.

1. Detection Reagents

The Oxidative Stress Detection Reagent (Green) is supplied lyophilized and should be reconstituted in 60 μ L anhydrous DMF to yield a 5 mM stock solution. Upon reconstitution, the stock solution should be stored at -20°C. The shelf life of the reconstituted reagent is about 1 week at -20°C. Gently mix before use.

The Superoxide Detection Reagent (Orange) is supplied lyophilized and should be reconstituted in 60 μ L anhydrous DMF to yield a 5 mM stock solution concentration. Upon reconstitution, the stock solution should be stored at -20°C. Gently mix before use.

2. Positive Control

The ROS Inducer (Pyocyanin) is supplied lyophilized and should be reconstituted in 20 μL anhydrous DMF to yield a 50 mM stock solution. For use, a final concentration of 200-500 μM is recommended. However, the optimal final concentration is cell-dependent and should be determined experimentally for each cell line being tested. ROS induction generally occurs within 20-30 minutes upon pyocyanin treatment and may decrease or disappear after that time. Plan accordingly.

3. Negative Control

The ROS Inhibitor (N-acetyl-L-cysteine) should be reconstituted in 123 μL of deionized water to yield a 0.5 M stock concentration. N-acetyl-cysteine is not readily soluble and may require vortexing. For use, a final concentration of 5 mM is recommended. However, the optimal final concentration is cell-dependent and should be determined experimentally for each cell line being tested.

Endogenous fluorescence of untreated cells should be determined in advance or per assay.

4. 1X Wash Buffer

Prepare 1X Wash Buffer by dissolving the contents of the pack in 1 liter of deionized water. When not in use, the 1X Wash Buffer should be stored refrigerated. Warm it to room temperature before use.

5. ROS/Superoxide Detection Solution (2X)

Prepare the ROS/Superoxide Detection Mix as follows: To every 10 mL of 1X Wash Buffer (see step 4) or culture medium (without phenol red), add 4 μL Oxidative Stress Detection Reagent (Green) and 4 μL Superoxide Detection Reagent (Orange). Gently mix to make 1:2500 diluted staining solution.

NOTE: For different cell lines, the staining efficiency might be different. 1:500 to 1:2500 is the suggested range for testing.

NOTE: Depending on the experiments, dyes can be used separately according to a provided protocol.

NOTE: To perform the compensation correction, single stained samples should be used. Make sure single component solutions (Green and Orange) are available.

B. CELL PREPARATIONS

Cells should be maintained via standard tissue culture practices. Always make sure that cells are healthy and in the log phase of growth before using them for the experiment.

C. FLUORESCENCE/CONFOCAL MICROSCOPY (ADHERENT CELLS)

1. The day before the experiment, seed the cells directly onto glass slides or polystyrene tissue culture plates to ensure ~ 50-70% confluency on the day of the experiment. It is important to have cells for both positive and negative controls.

IMPORTANT: Cells should be healthy and not overcrowded since results of the experiments will depend significantly on the cells' condition.

2. On the day of the experiments, change media to fresh media and label the wells as needed.
3. Negative control: add ROS inhibitor (N-acetyl-L-cysteine) to pre-treat the cells for at least 30 min prior to induction (see step A-3, page 5).
4. Induction: Load the same volume of the **ROS/Superoxide Detection Solution (2X)** (see step A-5, page 5) with addition of either the vehicle, experimental agent, ROS inducer (pyocyanin) at a desirable working concentration. Fully cover the cell monolayer. Incubate under normal tissue culture conditions for 30 min to 1 hour.

IMPORTANT: If the vehicle, experimental agent and ROS inducer (pyocyanin) will be added to the **ROS/Superoxide Detection Solution (2X)**, their concentration needs to be 2X. For the negative control, 1X NAC needs to be compensated in the solution to maintain its final concentration at the same level.

5. Carefully remove the ROS/Superoxide Detection Mix from the glass slides by gently tapping them against layers of paper towel, or from tissue culture plates.
6. Carefully wash cells twice with 1X Wash Buffer in a volume sufficient to cover the cell monolayer.
7. Add a few drops of 1X Wash Buffer on the top of the cells and immediately overlay the cells with a cover slip and observe them under a fluorescence/confocal microscope using standard excitation/emission filter sets. Oxidative stress detection requires a filter set compatible with Fluorescein (Ex/Em: 490/525 nm). Make sure prepared samples are protected from drying. Dried out cells may present different fluorescence patterns.

Recommended filter sets:

- Oxidative stress detection requires a filter set compatible with Fluorescein (Ex/Em: 490/525 nm).
- Superoxide detection requires a filter set compatible with Rhodamine (Ex/Em: 550/620 nm).

***Note:** Different exposure times may be required for optimal detection of the two dyes used in the kit.*

D. FLUORESCENCE/CONFOCAL MICROSCOPY (SUSPENSION CELLS)

1. Cells should be cultured to a density not to exceed 1×10^6 cells/ml. Make sure that cells are in the log phase of growth before starting an experiment.

***IMPORTANT:** Cells should be healthy and not overcrowded since results of the experiments will depend significantly on the cells' overall condition. A sufficient volume of cells should be centrifuged at $400 \times g$ for 5 minutes, yielding a working cell count of 1×10^5 cells/sample.*

2. Centrifuge cells for $400 \times g$ for 5 minutes and remove the supernatant.
3. Resuspend the cells in fresh media and aliquot 1×10^5 cells/sample at $1-5 \times 10^5$ cells/mL.
4. Negative control: add ROS inhibitor (N-acetyl-L-cysteine) to pre-treat the cells for at least 30 min prior to induction (see step A-3, page 5).

5. Induction: Load the same volume of the **ROS/Superoxide Detection Solution (2X)** (see step A-5, page 5) with addition of either the vehicle, experimental agent, ROS inducer (pyocyanin) at a desirable working concentration. Incubate under normal tissue culture conditions for 30 min to 1 hour with periodic shaking.

IMPORTANT: *If the vehicle, experimental agent and ROS inducer (pyocyanin) will be added to the **ROS/Superoxide Detection Solution (2X)**, their concentration needs to be 2X. For the negative control, 1X NAC needs to be compensated in the solution to maintain its final concentration at the same level.*

6. Centrifuge the cells at 400 x g for 5 minutes to remove the ROS Detection Solution.
7. Resuspend the cells in 5 mL of 1X Wash Buffer, centrifuge them at 400 x g for 5 minutes and remove the supernatant.
8. Resuspend the cells in 100 μ L of 1X Wash Buffer and apply a 20 μ L aliquot of the cell suspension, sufficient for 2×10^4 cells, onto a microscope slide. Immediately overlay the cells with a cover slip and analyze via fluorescence microscopy. Oxidative stress detection requires a filter set compatible with Fluorescein (Ex/Em: 490/525 nm). Make sure that prepared samples are protected from drying. Dried out cells may present different fluorescence patterns.

Recommended filter sets:

- Oxidative stress detection requires a filter set compatible with Fluorescein (Ex/Em: 490/525 nm).
- Superoxide detection requires a filter set compatible with Rhodamine (Ex/Em: 550/620 nm).

Note: *Different exposure times may be required for optimal detection of the two dyes used in the kit.*

E. FLOW CYTOMETRY (ADHERENT CELLS)

1. The day before the experiment, seed the cells on appropriate tissue culture plates to ensure ~ 50-70% confluency on the day of the experiment.

IMPORTANT: *Cells should be healthy and not overcrowded since results of the experiments will depend significantly on the cells' condition.*

2. Detach cells from the tissue culture plates using any appropriate method, collect cells in 5 mL round-bottom polystyrene tubes and wash them with 1X Wash Buffer. Centrifuge the cell suspension for 5 min. at 400 x g at room temperature. Discard the supernatant.
3. Simultaneously treat the cells with an experimental test agent (or control) and load the cells with the **ROS/Superoxide Detection Solution** (see step A-5, page 5).

i. Treatment conditions and controls

- a) Positive control samples should be established by treatment with ROS Inducer (Pyocyanin) (see step A-2, page 5)
- b) Negative control samples should be established by treatment with the ROS Inhibitor (N-acetyl-L-cysteine) (see step A-3, page 5)

NOTE: Cells should be pre-treated with the ROS Inhibitor at least 30 minutes prior to induction.

- c) Untreated samples should use the vehicle
- d) Experimental samples

ii. Loading the cells with dye

Resuspend cell pellet in 500 μ L of **ROS/Superoxide Detection Solution**, containing the treatment. Stain cells for 30 min. at 37°C in the dark. No washing is required prior to the analysis of the samples by flow cytometry.

IMPORTANT: Compensation correction will be needed to avoid overlap between green and orange fluorescent signals (see Appendix C).

Recommended controls for flow cytometry:

- Unstained untreated cells
- Positive control cells (pyocyanin-treated) stained with Oxidative Stress Detection Reagent (Green) only ("Green" cells)
- Positive control cells (pyocyanin-treated) stained with Superoxide Detection Reagent (Orange) only ("Orange" cells)

F. FLOW CYTOMETRY (SUSPENSION CELLS)

1. Cells should be cultured to a density not to exceed 1×10^6 cells/mL. Make sure that cells are in the log phase of growth before starting an experiment.

IMPORTANT: Cells should be healthy and not overcrowded since results of the experiments will depend significantly on the cells' overall condition. A sufficient volume of cells should be centrifuged at $400 \times g$ for 5 minutes, yielding a working cell count of $1-5 \times 10^5$ cells/sample.

2. Centrifuge cells for $400 \times g$ for 5 minutes and remove the supernatant.
3. Resuspend cells in fresh media at a concentration of $1-5 \times 10^5$ /mL. Aliquot 0.5-1 mL per sample into flow tubes.
4. Negative control: add ROS inhibitor (N-acetyl-L-cysteine) to pre-treat the cells for at least 30 min prior to induction (see step A-3, page 5).
5. Induction: Load the same volume of the **ROS/Superoxide Detection Solution (2X)** (see step A-5, page 5) with addition of either the vehicle, experimental agent, ROS inducer (pyocyanin) at a desirable working concentration. Incubate under normal tissue culture conditions for 30 min to 1 hour under dark with periodic shaking. No washing is required prior to the analysis of the samples by flow cytometry.

IMPORTANT: If the vehicle, experimental agent and ROS inducer (pyocyanin) will be added to the **ROS/Superoxide Detection Solution (2X)**, their concentration need to be 2X. For the negative control, 1X NAC needs to be compensated in the solution to maintain its final concentration at the same level.

IMPORTANT: Compensation correction will be needed to avoid overlap between green and orange fluorescent signals (see Appendix C).

Recommended controls for flow cytometry:

- Unstained untreated cells
- Positive control cells (pyocyanin-treated) stained with Oxidative Stress Detection Reagent (Green) only ("Green" cells)

- Positive control cells (pyocyanin-treated) stained with Superoxide Detection Reagent (Orange) only ("Orange" cells)

G. FLUORESCENCE MICROPLATE ASSAY (ADHERENT CELLS)

1. The day before the experiment, seed the cells in 96-well black wall/clear bottom plates at a density of $1-2 \times 10^4$ cells per well to ensure ~ 70-80% confluency on the day of the experiment. Leave several wells empty for the background fluorescence control measurements.

***Important:** Cells should be healthy and not overcrowded since results of the experiments will depend significantly on the cells' condition.*

2. Remove supernatant from cells.
3. Carefully wash cells with 1X Wash Buffer.
4. Simultaneously treat the cells with an experimental test agent (or controls) and load the cells with the **ROS/Superoxide Detection Solution** (see step A-5, page 5).

i. Treatment conditions and controls

For the highest quality data use at least 6 – 8 replicates for each condition tested.

- a) Positive control samples should be established by treatment with ROS Inducer (Pyocyanin) (see step A-2, page 5)
- b) Negative control samples should be established by treatment with the ROS Inhibitor (N-acetyl-L-cysteine) (see step A-3, page 5)
***NOTE:** Cells should be pre-treated with the ROS Inhibitor at least 30 minutes prior to induction.*
- c) Untreated samples should use the vehicle
- d) Experimental samples

ii. Loading the cells with dye

Add 100 µL/well of **ROS/Superoxide Detection Solution**, containing the treatment. Stain cells for 60 min. at 37°C in the dark.

5. Read the plates (bottom reading), without removing the detection mix, using a fluorescence microplate reader and standard fluorescein (Ex=488 nm, Em=520 nm) and rhodamine (Ex=550 nm, Em=610 nm) filter sets.
6. If required, Z' factor may be calculated for each detection profile using the following formula:(5)

$$Z' = 1 - [(3 * SD_{\text{sample}} + 3 * SD_{\text{control}}) / (|Mean_{\text{sample}} - Mean_{\text{control}}|)]$$

H. FLUORESCENCE MICROPLATE ASSAY (SUSPENSION CELLS)

NOTE: Perform all steps requiring centrifugation at 400 x g for 5 minutes using a centrifuge with swinging buckets.

1. Collect the cells by centrifugation at 400 x g for 5 min, resuspend them in the appropriate cell culture medium (without phenol red) at a density of 0.5 x 10⁶ - 1.0 x 10⁶ cells/mL, count and aliquot 100 µL of the suspension into wells of 96-well black wall plates. Leave several wells empty for the background fluorescence control measurements.

Important: Cells should be healthy and not overcrowded since results of the experiments will depend significantly on the cells' condition.

2. Remove the medium by centrifugation of the plate.
3. Wash cells by centrifugation with 1X Wash Buffer.
4. Simultaneously treat the cells with an experimental test agent (or controls) and load the cells with the **ROS/Superoxide Detection Solution** (see step A-5, page 5).

iii. Treatment conditions and controls

For the highest quality data use at least 6 – 8 replicates for each condition tested.

- a) Positive control samples should be established by treatment with ROS Inducer (Pyocyanin) (see step A-2, page 5)

- b) Negative control samples should be established by treatment with the ROS Inhibitor (N-acetyl-L-cysteine) (see step A-3, page 5)

NOTE: *Cells should be pre-treated with the ROS Inhibitor at least 30 minutes prior to induction.*

- c) Untreated samples should use the vehicle
- d) Experimental samples

iv. Loading the cells with dye

Add 100 µL/well of **ROS/Superoxide Detection Solution**, containing the treatment. Stain cells for 60 min. at 37°C in the dark.

- 5. Read the plates (bottom reading), without removing the detection mix, using a fluorescence microplate reader and standard fluorescein (Ex=488 nm, Em=520 nm) and rhodamine (Ex=550 nm, Em=610 nm) filter sets
- 6. If required, Z' factor may be calculated for each detection profile using the following formula:(5)

$$Z' = 1 - [(3*SD_{\text{sample}}+3*SD_{\text{control}})/(|\text{Mean}_{\text{sample}} - \text{Mean}_{\text{control}}|)]$$

VI. APPENDICES

A. FILTER SET SELECTION

For fluorescence microscopy, careful consideration must be paid to the selection of filters. Dichroic filters should be selected in which the “cut-off” frequency is optimally mid-way between the two emission bands that are desired (one reflected, the other transmitted). However, it is important to realize that dichroic filters have a somewhat limited reflectance range, *i.e.*, a 600 nm short-pass dichroic filter may actually reflect light <500 nm. When selecting filters, it is critical to discuss with the filter or microscope manufacturer exactly what wavelength specifications are required for both the transmitted and the reflected light.

In addition, filters should be obtained that have the highest possible transmission efficiency (typically requiring anti-reflection coating). Each optic that an emission beam must traverse will remove some fraction of the desired light. The difference between 80% transmission and 95% transmission for each detector may result in a greater than three-fold difference in the amount of light available to the detector.

B. SETTING UP OPTIMAL EXPOSURE TIME FOR DETECTION OF THE DYE

Optimal exposure times should be established experimentally for each dye used in the experiment. Both negative and positive controls should be utilized. Start with the negative control (untreated stained cells) and set up the exposure time so the fluorescent background is negligible. Then switch to a positive control (pyocyanin-treated cells) and adjust the exposure time to record a bright fluorescent image. Avoid saturation of the signal (very bright spots on the image). If saturation of the signal occurs, decrease the exposure time. It is recommended to acquire 5-6 single color images for each sample.

C. COMPENSATION CORRECTION FOR FLOW CYTOMETRY

1. Signals produced by peroxides, peroxynitrite and hydroxyl radicals will be detected in the FL1 channel. Superoxide production will be detected in the FL2 channel. To avoid overlap between green and orange fluorescent signals the following compensation procedure should be performed.
2. Run the unstained uninduced sample first. Generate a FSC versus SSC dot plot and gate out cell debris.
3. Generate a log FL1 (X-axis) versus a log FL2 (Y-axis) dot plot. Adjust PMT voltages for both channels so the signals from unstained cells should fall within the first log decade scale of FL1 and FL2 axes.
4. Run single stained “Green” positive control and adjust FL2-%FL1 compensation until the orange fluorescence signal will fall into the first decade of the log FL2 scale.
5. Repeat compensation procedure with the “Orange” single stained positive control and adjust FL1-%FL2 compensation until the green fluorescence signal will fall into the first decade of the log FL1 scale.

NOTE: It is important to use the brightest positive single stained samples for proper compensation correction that allows distinguishing between negative and slightly positive (dim) cells.

D. ANTICIPATED RESULTS (FLUORESCENCE MICROSCOPY)

1. The Superoxide Detection Reagent (Orange) yields an evenly distributed, bright orange nuclear staining pattern in induced cells. Note the structural change in positively treated cells versus control untreated cells (diffuse, dim cytoplasmic structural pattern observed in the control cells is replaced with uniform cytoplasmic staining and bright nuclear staining in superoxide-positive cells).
2. Increased levels of oxidative stress give a uniform green cytoplasmic staining in the presence of the Oxidative Stress Detection Reagent (Green).
3. ROS positive control samples, induced with ROS Inducer (Pyocyanin), exhibit a bright orange fluorescence in the nucleus as well as a bright green fluorescence in the cytoplasm.
4. Cells pretreated with the ROS Inhibitor (N-acetyl-L-cysteine) should not demonstrate any green or orange fluorescence upon induction.
5. Untreated samples should present only low autofluorescent background signal in any channel.

E. FLOW CYTOMETRY DATA ANALYSIS AND ANTICIPATED RESULTS

1. It is critical that positive (pyocyanin-induced) and control (untreated) samples be included in every experiment for every cell type. Negative (NAC-pretreated) sample is optional but very helpful. In preliminary experiments, it is important to establish appropriate doses of inducers and inhibitors for each cell type used.
2. Cell debris should be gated out using FSC versus SSC dot plot.
3. Generate a log FL1 (X-axis) versus a log FL2 (Y-axis) dot plot and add quadrants to it. Adjust quadrants so the majority of control cells (80-90%) will fall into lower left quadrant. Keep the same quadrant gate throughout the assay.

NOTE: Remember that different cell types demonstrate different redox profiles therefore the number of the cells in the lower left quadrant may vary significantly between the cell lines.

4. Cells with increased production of superoxide demonstrate bright orange fluorescence and will be detected using the FL2 channel. Such cells will appear in the two upper quadrants of a log FL1 (X-axis) versus a log FL2 (Y-axis) dot plot.
5. Cells with increased levels of oxidative stress demonstrate a bright green staining in the presence of the Oxidative Stress Detection Reagent and can be registered in FL1 channel. Such cells will appear in the upper and lower right quadrants of a log FL1 (X-axis) versus a log FL2 (Y-axis) dot plot.
6. ROS positive control samples, induced with ROS Inducer (Pyocyanin), exhibit both bright orange and green fluorescence and appear to be positive in FL1 and FL2 channels. The increase of the cell population in the upper left, upper right and lower right quadrants will be registered.
7. Cells pretreated with the ROS Inhibitor (N-acetyl-L-cysteine) should not demonstrate significant green or orange fluorescence upon induction.

8. Control (untreated) samples should present only low auto-fluorescent background signal in any channel thus falling into the lower left quadrant on an FL1 versus FL2 dot plot.
9. Results of the experiments can be presented as percentage of the cells with increased ROS production or as increase in the mean fluorescence of the induced samples versus control.

F. FLUORESCENCE MICROPLATE ASSAY DATA ANALYSIS AND ANTICIPATED RESULTS

1. It is critical that positive (pyocyanin-induced) and control (untreated) samples be included in every experiment for every cell type. Negative (NAC-pretreated) sample is optional, but very helpful. In preliminary experiments, it is important to establish appropriate doses of inducers and inhibitors for each cell type used.
2. Cells with increased production of superoxide demonstrate bright orange fluorescence in the presence of the Superoxide Detection Reagent and the signal will be detected using standard rhodamine filter settings.
3. Cells with increased levels of oxidative stress demonstrate a bright green staining in the presence of the Oxidative Stress Detection Reagent and the signal will be detected using standard fluorescein filter setting.
4. ROS positive control samples, induced with ROS Inducer (Pyocyanin), exhibit both bright orange and green fluorescence and can be detected in both green and orange channels. Cells pretreated with the ROS Inhibitor (N-acetyl-L-cysteine) should not demonstrate significant green or orange fluorescence upon induction.
5. Control (untreated) samples should present only low autofluorescence signal in both channels.

NOTE: Remember that different cell types demonstrate different redox profiles. Therefore, the autofluorescence signal may vary significantly in both channels.

6. Results of the experiments should be normalized using the background readings from empty wells (see step H1) and can be presented as ratios of the mean fluorescence of the induced samples versus control for each channel.

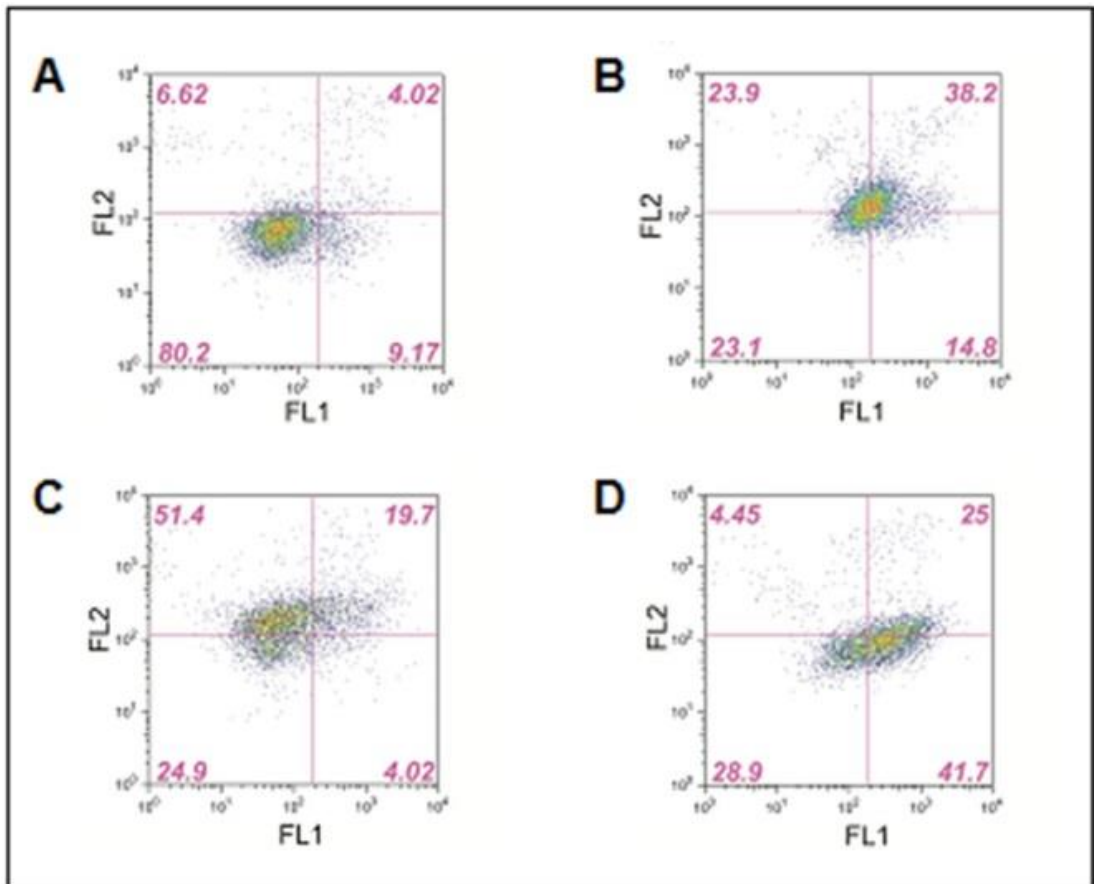


Figure 1. Jurkat cells were induced with 100 μ M pyocyanin (general ROS inducer, panel B), 200 μ M antimycin A (superoxide inducer, panel C) or 1 μ M of t-butyl-hydroperoxide (peroxide inducer, panel D), stained with two color ROS Detection Kit and analyzed using flow cytometry. Untreated cells (panel A) were used as a control. Cell debris were ungated and compensation was performed using single stained pyocyanin-treated samples. Red numbers reflect the percentage of the cells in each quadrant.

VII. REFERENCES

1. Tarpey, M. and Fridovich, I. *Circ Res.* 89 (2001), 224-236.
2. Batandier, C., et al. *J Cell Mol Med.* 6 (2002), 175-187.
3. Gomes, A., et al. *J Biochem Biophys Meth.* 65 (2005), 45-80.
4. Wardman, P. *Free Rad Biol Med.* 43 (2007), 995-1022.
5. Zhang, J.H., et al. *J Bio Screening* 4 (1999).

VIII. TROUBLESHOOTING GUIDE

Problem	Potential Cause	Suggestion
Low or no fluorescent signal in positive control	Dead or stressed (overcrowded) cells	Prepare fresh cell culture for the experiments. Make sure that the cells are in the log growth phase.
	Band pass filters are too narrow or not optimal for fluorescent probes (fluorescence microscopy)	Multiple band pass filters sets provide less light than single band pass ones. Check Methods and Procedures section of this manual and Appendix A for recommendations.
	Insufficient fluorescent dye concentration	Follow the procedures provided in this manual.
	Insufficient inducer concentration	Determine an appropriate concentration of inducer for the cell line(s) used in the study.
	Species of interest may react with each other, thus attenuating the expected signal.	Check signaling pathways and all the components present in the cellular environment.
	Inappropriate time point of the detection	Make sure that time of detection is optimized and the samples are prepared immediately. Green signal may quench if concentration of product becomes too high (due to long exposure to the inducer). Otherwise, oxidized product may eventually leak out of the cells when left for a prolonged period.
	Overcompensation of the signal (flow cytometry)	Change the values of compensation correction using single stained positive samples. Follow recommendation in the Appendix.

Problem	Potential Cause	Suggestion
Low or no fluorescent signal in positive control	Cell density is too low in microplate well.	Check the cell count to confirm proper cell density. For suspension cells, careful removal of supernatant after washing steps is critical, as cells may be dislodged and washed away.
High fluorescent background	Stressed (overcrowded) cells	Prepare new cell culture for the experiment. Make sure that the cells are in the log growth phase.
	Band pass filters are too narrow or not optimal for fluorescent probes.	Use correct filter for each fluorophore. Check Methods and Procedures section of this manual and Appendix A for the recommendations.
	Wash step is necessary.	Follow the procedures provided in this manual, making optional wash steps mandatory.
	Inappropriate time point for detection	Make sure that time of detection is optimized and the samples are prepared immediately.
	Inappropriate cell conditions	Make sure that you have viable cells at the beginning of the experiment, and that the inducer treatment does not kill the cells during the time frame of the experiment.
No decrease in the fluorescent signal after using a specific inhibitor	Inappropriate inhibitor concentration (too low or too high)	Very low doses of inhibitor may not affect ROS production by inducer. Alternatively, very high doses of the inhibitors may cause oxidative stress itself and generate fluorescent signal. Optimize the concentration of the inhibitor and pretreatment time for each particular cell line.

Problem	Potential Cause	Suggestion
No decrease in the fluorescent signal after using a specific inhibitor	Inappropriate time point for detection	When cells are kept too long with the inhibitors or at very high inducer concentrations, after a certain time, the inhibitor becomes insufficient. Make sure that time of detection is optimized.
	Inappropriate filter set on the microscope	Use correct filter for excitation and emission. Check Methods and Procedures section of this manual and Appendix A for the recommendations. Minimal spectral overlap should occur with the selected set of filters.



Product Manual

NOTES



Product Manual

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