



Product Manual

ROS-ID[®] ROS/RNS Detection Kit

for fluorescence microscopy

200 assays

Instruction Manual

Catalog Number: **ENZ-51001-200**



Product Manual

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.

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.....	7
Fluorescence/Confocal Microscopy (Adherent cells).....	7
Fluorescence/Confocal Microscopy (Suspension cells).....	9
Appendices.....	11
Filter Set Selection.....	11
Setting up Optimal Exposure Time, Detection of the Dye.....	11
Anticipated Results.....	11
References.....	12
Troubleshooting Guide.....	13
Contact Information.....	18



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



Please contact Enzo Life Sciences Technical Support if necessary.

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[®] ROS/RNS Detection Kit enables detection of comparative levels of ROS/RNS in cells and also distinguishes between different reactive species in live cells. Through the combination of three specific fluorescent probes, the kit provides a simple and specific assay for the real-time measurement of free nitric oxide (NO) and by extension nitric oxide synthase (NOS) activity, as well as global levels of reactive oxygen species (ROS), and specifically superoxide in living cells.

This kit is designed to directly monitor real time reactive oxygen and/or nitrogen species (ROS/RNS) production in live cells using fluorescence microscopy. The kit includes three fluorescent dye reagents as major components: NO Detection Reagent (red fluorescent), Oxidative Stress Detection Reagent (green fluorescent) for total ROS detection reagent and Superoxide Detection Reagent (orange fluorescent). The non-fluorescent, cell-permeable NO detection dye reacts with NO in the presence of O₂ with high specificity, sensitivity and accuracy, yielding a water-insoluble red fluorescent product. The non-fluorescent, cell-permeable total ROS detection dye reacts directly with a wide range of reactive species, such as hydrogen peroxide, peroxynitrite and hydroxyl radicals, yielding a green fluorescent product indicative of cellular production of different ROS/RNS types. The superoxide detection dye is a cell-permeable probe that reacts specifically with superoxide, 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 three dyes can be visualized using a wide-field fluorescence microscope equipped with standard green (490/525 nm), orange (550/620 nm), and red (650/670 nm) fluorescent cubes.

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 assays using live cells (adherent or in suspension)

Reagent	Quantity
NO Detection Reagent(Red)	60 µL
Oxidative Stress Detection Reagent (Green)	300 nmoles
Superoxide Detection Reagent (Orange)	300 nmoles
NO Inducer (L-Arginine)	100 µL
ROS Inducer (Pyocyanin)	1 µmole
NO Scavenger (c-PTIO)	400 nmoles
ROS Inhibitor (N-acetyl-L-cysteine)	2 x 10 mg
10X Wash Buffer	15 mL

III. OTHER MATERIALS NEEDED

- CO₂ incubator (37°C)
- Standard fluorescence microscope
- Calibrated, adjustable precision pipettors, preferably with disposable plastic tips
- Tubes appropriate for holding cells during induction of ROS/RNS (for suspension cells only)
- 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

- This product is for research use only and is not intended for diagnostic purposes.
- The NO Detection Reagent (Red) contains DMF 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.



Protect
from light

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

1. Detection Reagents

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

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

2. Positive Control

- The NO Inducer (L-Arginine) is supplied at a stock concentration of 100 mM in deionized water. A final concentration of 1 mM is recommended. However, the optimal final concentration is cell-dependent and should be determined experimentally for each cell line being tested. NO induction generally occurs within 15-20 minutes upon treatment with NO Inducer (L-Arginine) and may drastically decrease or disappear after that time. Plan accordingly.
- The ROS Inducer (Pyocyanin) is supplied lyophilized and should be reconstituted in 100 μ L anhydrous DMF to yield a 10 mM stock solution. For use, a final concentration of 200-500 μ M 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 NO Scavenger (c-PTIO) is supplied lyophilized and should be reconstituted in 100 μ L DMF to produce a stock solution of 4 mM. For use, a final concentration of 20-80 μ M is recommended. However, the optimal final concentration is cell-dependent and should be determined experimentally for each cell line being tested.

Note: *Adherent cells pretreated with c-PTIO may become weakly adherent and/or detach from the cell culture substratum.*

- The ROS Inhibitor (N-acetyl-L-cysteine) should be reconstituted in 123 μ L of deionized water to yield a 0.5 M stock solution. 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

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

5. ROS-ID[®] ROS/RNS 3-Plex Detection Mix

- Pre-dilute the reconstituted Oxidative Stress Detection Reagent (Green) stock 1:10 using pre-warmed tissue culture media.
- Pre-dilute the reconstituted Superoxide Detection Reagent (Orange) stock 1:10 using pre-warmed tissue culture media.

Prepare the ROS-ID[®] ROS/RNS 3-Plex Detection Mix by combining appropriate volumes of the pre-diluted Oxidative Stress Detection Reagent (step V.A.1, page 3), pre-diluted Superoxide Detection Reagent (step V.A.2, page 4) and NO Detection Reagent (Red) using the volumes specified in the following table.

Note: *The NO Detection Reagent (Red) is ready-to-use and does not require a pre-dilution step.*

Important: *Pre-warmed media should be used to avoid precipitation of the NO Detection Reagent (Red) in the mix.*

ROS-ID® ROS/RNS 3-Plex Detection Mix	
Reagent	Vol. Req'd per 1 mL Detection Mix
NO Detection Reagent(Red)	2.5 µL
Diluted (1:10) Oxidative Stress Detection Reagent (Green) from step V.A.1	2 µL
Diluted (1:10) Superoxide Detection Reagent (Orange) from step V.A.2	2 µL
Tissue Culture Medium	993.5 µL
Total Volume	1 mL

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

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.

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

Load the cells with the ROS-ID® ROS/RNS 3-Plex Detection Mix using a volume sufficient to cover the cell monolayer and incubate under normal tissue culture conditions for 2 hours.

Carefully remove the ROS-ID® ROS/RNS 3-Plex Detection Mix from the glass slides by gently tapping them against layers of paper towel, or from tissue culture plates. **Optional:** Cells may be washed with the 1X Wash Buffer.

2. Treat the cells with an experimental test agent. Separate positive control samples should be treated with the NO Inducer (L-Arginine) and the ROS Inducer (Pyocyanin). Negative control samples should be established by treatment with the ROS Inhibitor (N-acetyl-L-cysteine) or NO Scavenger (c-PTIO) correspondingly, for the same length of time as the inducer.

Note: *Cells should be treated with the NO Scavenger or ROS Inhibitor 30 minutes prior to induction.*

All treatments should be performed under normal tissue culture conditions. It is recommended to perform a pretreatment by adding the NO Scavenger or ROS Inhibitor to the aliquots of ROS-ID[®] ROS/RNS 3-Plex Detection Mix for the last 30 minutes of the reagent loading. Treatment with an experimental test agent or ROS/NO inducers included with the kit should be performed in the cell culture media without dyes.

3. Carefully wash cells twice with 1X Wash Buffer in a volume sufficient to cover the cell monolayer.

Note: *Adherent cells pre-treated with NO Scavenger (c-PTIO) may become loose and/or detach from the cell culture substratum.*

4. Overlay the cells with a cover slip and observe them under a fluorescence/confocal microscope using standard excitation/emission filter sets. 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).
- NO detection requires a filter set compatible with Cyanine 5 (650/670 nm).

Note: *Different exposure times may be required for optimal detection of the three 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.

Resuspend the cell pellet in 100 μ L of ROS-ID[®] ROS/RNS 3-Plex Detection Mix (see step V.A.5, page 5) and incubate under normal tissue culture conditions for 2 hours with periodic shaking.

Centrifuge the cells at $400 \times g$ for 5 minutes to remove the ROS-ID[®] ROS/RNS 3-Plex Detection Mix. Optional: Resuspend the cells in 0.5 mL 1X Wash Buffer, centrifuge them at $400 \times g$ for 5 minutes and remove the supernatant.

2. Treat the cells with an experimental test agent. Separate positive control samples should be treated with the NO Inducer (L-Arginine) and the ROS Inducer (Pyocyanin). Negative control samples should be established by treatment with the ROS Inhibitor (N-acetyl-L-cysteine) or NO Scavenger (c-PTIO) correspondingly, for the same length of time as the inducer.

Note: Cells should be treated with the NO Scavenger or ROS Inhibitor 30 minutes prior to induction.

All treatments should be performed under normal tissue culture conditions. It is recommended to perform a pretreatment by adding the NO Scavenger or ROS Inhibitor to the aliquots of ROS-ID[®] ROS/RNS 3-Plex Detection Mix for the last 30 minutes of the reagent loading. Treatment with an experimental test agent or ROS/NO inducers included with the kit should be performed in the cell culture media without dyes.

3. Centrifuge the cells at $400 \times g$ for 5 minutes.

4. Resuspend the cells in 0.5 mL of 1X Wash Buffer, centrifuge them at 400 x g for 5 minutes and remove the supernatant.
5. 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. Overlay the cells with a cover slip and analyze immediately via fluorescence microscopy. 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).
- NO detection requires a filter set compatible with Cyanine 5 (650/670 nm).

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

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. ANTICIPATED RESULTS

1. In the presence of NO, the NO Detection Reagent (Red) will demonstrate a red punctate cytoplasmic staining pattern.
2. 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).

3. Increased levels of oxidative stress give a uniform green cytoplasmic staining in the presence of the Oxidative Stress Detection Reagent.
4. Nitric oxide (NO) positive control samples, induced with NO Inducer (L-Arginine), should exhibit red fluorescence, with punctate cytoplasmic staining pattern. Keep in mind that fluorescence of the NO Detection Reagent (Red) is not very bright. Depending on the filters used, the spill-over of the red signal may be visible in the orange channel. However, in this case, the staining pattern will differ significantly from the superoxide generated signal (punctate cytoplasmic staining versus nuclear staining). Moreover, this staining should disappear upon pretreatment with NO Scavenger (c-PTIO).
5. 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.
6. Cells pretreated with the ROS Inhibitor (N-acetyl-L-cysteine) should not demonstrate any green or orange fluorescence upon induction.
7. Cells incubated with the NO Scavenger (c-PTIO), post-induced with NO Inducer (L-Arginine) should not demonstrate red fluorescence
8. Untreated samples should present only low auto fluorescent background signal in any channel.

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.

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.	Multiple band pass filters sets provide less light than single band pass ones. Use correct filter for each fluorophore. Check Methods and Procedures section of this manual and Appendix A for recommendations. Use of a wide band pass or a long pass filters is recommended (in particular, for the NO detection).
	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. Red signal may disappear over time because the NO oxidized fluorescent probe can undergo further redox changes in the cellular physiological environment. 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.

Problem	Potential Cause	Suggestion
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. Minimal spectral overlap should occur with the selected set of filters.
	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 of 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.
	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.



Product Manual

Problem	Potential Cause	Suggestion
	Inappropriate filter set on the microscope	Use correct filter for each fluorophore. 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

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