

ROS-ID[®] NO Detection kit

Enzo Life Sciences' ROS-ID[®] NO Detection kit is designed to directly monitor real time NO production in live cells by fluorescence microscopy using NO Detection Reagent (red fluorescent) as the major component. 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. Importantly, this dye is not reactive towards peroxynitrite that allows distinguishing between peroxynitrite and nitric oxide. 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 this dye can be visualized using a wide-field fluorescence microscope equipped with any red fluorescent cube. The combination of 650/670 nm is recommended when additional fluorescence signals (green or orange) would be detected simultaneously. The NO Detection Kit contains sufficient reagents for at least 200 microscopy assays using live cells (adherent or in suspension).

Citations: 24

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-51013-200	1Kit
---------------	------

Manuals, SDS & CofA

[View Online »](#)

- Directly monitor real time NO production in live cells by fluorescence microscopy
- High specificity, sensitivity and accuracy, yielding a water-insoluble red fluorescent product
- Distinguish between peroxynitrite and nitric oxide
- NO inducer and scavenger included

Handling & Storage

Handling	Protect from light. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Fluorescence microscopy, Fluorescent detection
Application Notes	This kit is designed to directly monitor real time NO production in live cells using fluorescence microscopy.
Contents	NO Detection Reagent (Red), 60 µl NO Inducer (L-Arginine), 100 µl NO Scavenger (c-PTIO), 400 nmoles 10X Wash Buffer, 15 ml
Quality Control	<p>A sample from each lot of ROS-ID[®] NO Detection kit is used to stain HeLa cells using the procedures described in the user manual. The stained cells are analyzed using a wide-field fluorescence microscope equipped with standard red (650/670 nm) fluorescent cubes.</p> <p>The following results are obtained: Nitric Oxide positive control samples induced with L-Arginine exhibit red fluorescence with a red punctuate cytoplasmic staining pattern. Cells incubated with the NO Scavenger (c-PTIO), post induced with L-Arginine don't demonstrate any red fluorescence.</p>
Quantity	200 fluorescence microscopy assays.
Technical Info / Product Notes	<p>The ROS-ID[®] NO detection kit is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.</p>



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)