

# MITO-ID<sup>®</sup> Red detection kit (GFP- CERTIFIED<sup>®</sup>)

## Photostable, non-toxic and selective mitochondrial dye that stains regardless of

**Membrane potential**  
MITO-ID<sup>®</sup> Red Detection Kit (GFP-CERTIFIED<sup>®</sup>) contains a proprietary membrane-permeable mitochondria-selective dye suitable for use with live-, detergent-permeabilized- and even aldehyde-fixed-cells. Unlike conventional dyes, such as DiOC6(3), JC-1, rhodamine 123 and tetramethylrhodamine ethyl ester, MITO-ID<sup>®</sup> Red dye highlights mitochondria regardless of their energetic state. The dye is compatible with most fluorescence detection systems, including conventional and confocal fluorescence microscopes, as well as, High Content Screening (HCS) platforms. The kit is useful for assessing mitochondrial morphology changes, estimating mitochondrial mass and co-localizing GFP-tagged proteins to the mitochondrial compartment. This kit is specifically designed for use with GFP expressing cell lines, as well as cells expressing blue, cyan or yellow fluorescent proteins (BFPs, CFPs, YFPs). Additionally, the kit is suitable for use with live or post-fixed cells in conjunction with fluorescent probes, such as labeled antibodies, or other fluorescent conjugates displaying similar spectral properties such as fluorescein and coumarin. A nuclear counterstain (Hoechst 33342) is provided to highlight this organelle as well. Wavelength maxima: MITO-ID<sup>®</sup> Red  $\lambda_{ex}$  558 nm,  $\lambda_{em}$  690 nm; Hoechst 33342  $\lambda_{ex}$  350 nm,  $\lambda_{em}$  461 nm.

Citations: 38

[View Online »](#)

## Ordering Information

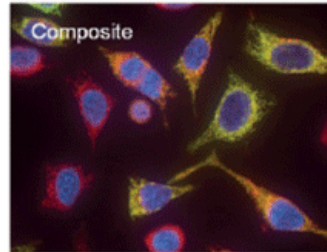
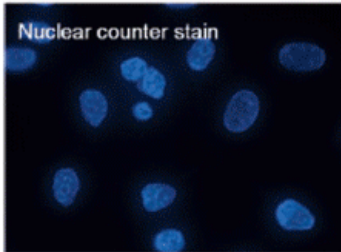
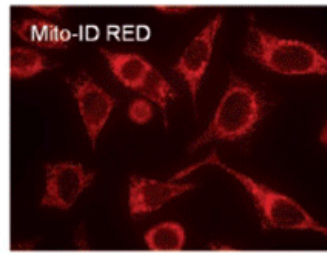
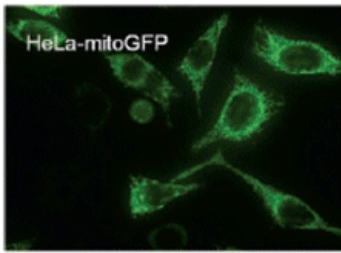
[Order Online »](#)

ENZ-51007-0100	100 tests
ENZ-51007-500	500 tests

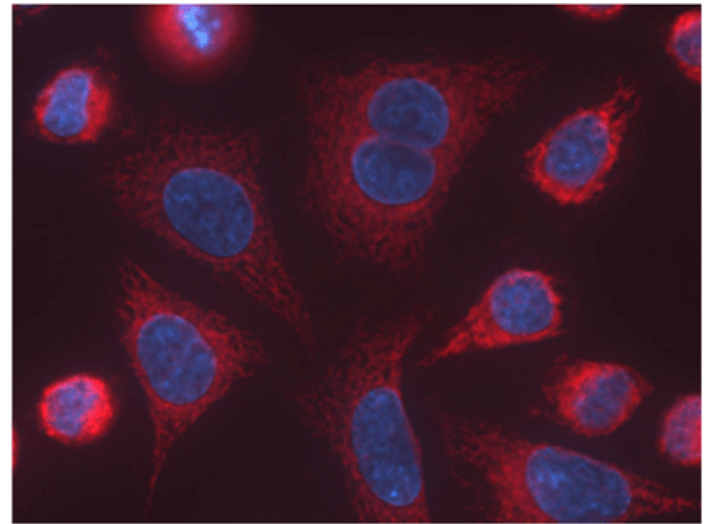
## Manuals, SDS & CofA

[View Online »](#)

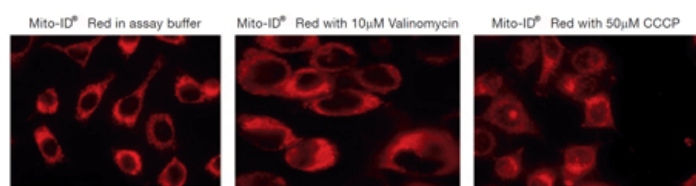
- Mitochondria-selective dye stains live, detergent permeabilized and aldehyde fixed cells
- Long wavelength red emission is easily multiplexed with common fluorescent dyes
- Highly resistant to photobleaching and concentration quenching, for strong, consistent fluorescent signal
- Highlights mitochondria regardless of the organelle's membrane potential status
- Stringently manufactured, to control and eliminate non-specific assay artifacts



HeLa-TurboGreen-mitochondria cells (HeLa-mitoGFP, MarinPharm GmbH, Luckenwalde, Germany) stained with MITO-ID<sup>®</sup> Red and Hoechst 33342 (blue) dyes. MITO-ID<sup>®</sup> Red co-localizes with the EGFP-cytochrome C oxidase signal (yellow signal), demonstrating selectivity for mitochondria. Note that mitochondria in cells no longer expressing the GFP-tagged protein appear red in the composite image.



Composite fluorescence microscopy images of HeLa cells (40X objective lens). Cells were stained with MITO-ID<sup>®</sup> Red dye for 15 minutes. Nuclei were counter-stained with Hoechst 33342 dye.



The MITO-ID<sup>®</sup> Red 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 perturbing drugs, such as valinomycin.



MITO-ID<sup>®</sup> Red detection kit (GFP-CERTIFIED<sup>®</sup>)

## Handling & Storage

Use/Stability	With proper storage, the kit components are stable up to the date noted on the product label. Store kit at -20°C in a non-frost free freezer, or -80°C for longer term storage.
Handling	Protect from light. Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	Fluorescence microscopy, Fluorescent detection
Application Notes	For use with GFP-expressing cell lines, as well as cells expressing blue, cyan or yellow fluorescent proteins.
Contents	MITO-ID <sup>®</sup> Red Detection Reagent Hoechst 33342 Nuclear Stain 10X Assay Buffer
Quality Control	<ol style="list-style-type: none"><li>1. Absorption peak of MITO-ID<sup>®</sup> Red Detection Reagent: <math>\lambda_{\text{max}} = 550 \pm 7 \text{ nm}</math></li><li>2. % purity of MITO-ID<sup>®</sup> Red Detection Reagent by HPLC: <math>\geq 93\%</math></li><li>3. A sample from each lot of GFP-CERTIFIED<sup>®</sup> MITO-ID<sup>®</sup> Red Mitochondrial Detection Kit is used to stain HeLa cells, expressing GFP-cytochrome C oxidase, using the procedures described in the user manual. The selectivity of the MITO-ID<sup>®</sup> Red dye is evident as shown by the co-localization with a GFP-cytochrome C oxidase fusion expressed in the HeLa cells.</li></ol>
Quantity	For ENZ-51007-500, 500 assays For ENZ-51007-0100, 100 assays

## Technical Info / Product Notes

### **Application Note:**

[Towards Understanding the Molecular Basis of Parkinson's Disease: Cell-based Model of Mitophagy and Aggresome Accumulation](#)



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)