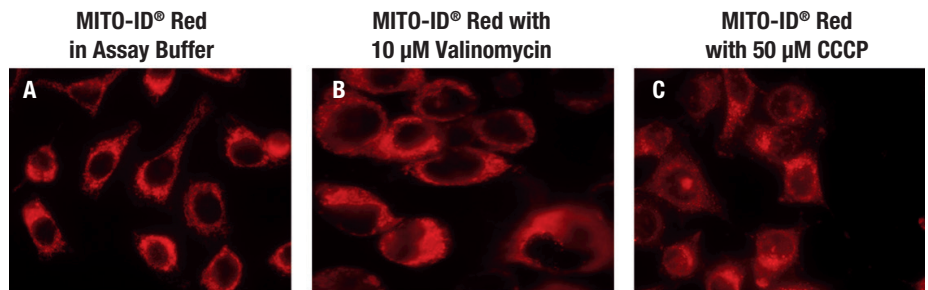


MITO-ID® Red & Green Detection Kits (ENZ-51007, ENZ-51022)

Photostable, Non-Toxic and Selective Mitochondrial Dye that Stains Regardless of Membrane Potential

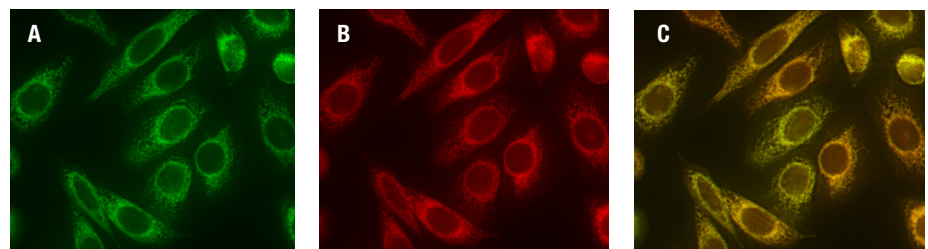
MITO-ID® Red (GFP-CERTIFIED®) and MITO-ID® Green Detection Kits contain proprietary membrane-permeable, mitochondria-selective dyes suitable for use with live, detergent-permeabilized and aldehyde-fixed cells. The kits are useful in assessing mitochondrial morphology changes, estimating mitochondrial mass, and co-localizing RFP-tagged proteins to the mitochondrial compartment. A blue nuclear counterstain, Hoechst 33342, is provided to highlight the nucleus.

- Potent mitochondria-selective dye suitable for live, permeabilized, and fixed cell staining
- Easily multiplexed with common fluorescent dyes and proteins
- Highly resistant to photobleaching, concentration quenching, and photo-conversion
- Highlights mitochondria regardless of the organelle's membrane potential status
- Stringently manufactured to control and eliminate non-specific assay artifacts



The MITO-ID® 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.

HeLa Cells Stained with MITO-ID® Green and Red Dyes



HeLa cells were stained with both MITO-ID® Green (A) and MITO-ID® Red (B) dyes. MITO-ID® Green dye co-localizes with the MITO-ID® Red dye, shown as yellow signal (C), demonstrating selectivity for mitochondria. Previously, MITO-ID® Red dye has been shown to co-localize with mitochondria expressing GFP.

RELATED PRODUCTS

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
DiOC6(3) iodide (Ultra Pure)	ENZ-52303	LYSO-ID® Red Detection Kit (GFP-CERTIFIED®)	ENZ-51005
ER-ID® Green Assay Kit	ENZ-51025	MITO-ID® Membrane Potential Cytotoxicity Kit	ENZ-51019
JC-1 (Ultra Pure)	ENZ-52304	MITO-ID® Membrane Potential Detection Kit	ENZ-51018
JC-10 (Ultra Pure)	ENZ-52305	Rhodamine 123 (Ultra Pure)	ENZ-52307
LYSO-ID® Green Detection Kit	ENZ-51034	TMRE (Ultra Pure)	ENZ-52309

Global Headquarters
ENZO LIFE SCIENCES, INC.
 Ph: 800.942.0430
 info-usa@enzolifesciences.com

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

Belgium, The Netherlands & Luxembourg
& Luxembourg
 Ph: +32 3 466 0420
 info-be@enzolifesciences.com

France
 Ph: +33 472 440 655
 info-fr@enzolifesciences.com

Germany
 Ph: +49 7621 5500 526
 info-de@enzolifesciences.com

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