

ORGANELLE-ID-RGB[®] reagent I

Mixture of fluorescent dyes for detection of lysosomes, mitochondria and nucleic acids

The ORGANELLE-ID-RGB[®] reagent I is a mixture of cell-permeable red fluorescent lysosomal dye, green fluorescent mitochondrial dye and blue fluorescent nucleic acid dye. The staining pattern arising from the combination of these three dyes permits visualization of the target organelles by fluorescence/confocal microscopy. The reagent, supplied as a 500X solution, is sufficient for 1000 microscopy assays. The single-tube format makes this multi-organelle stain reagent easy to use.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-53007-C200	200µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

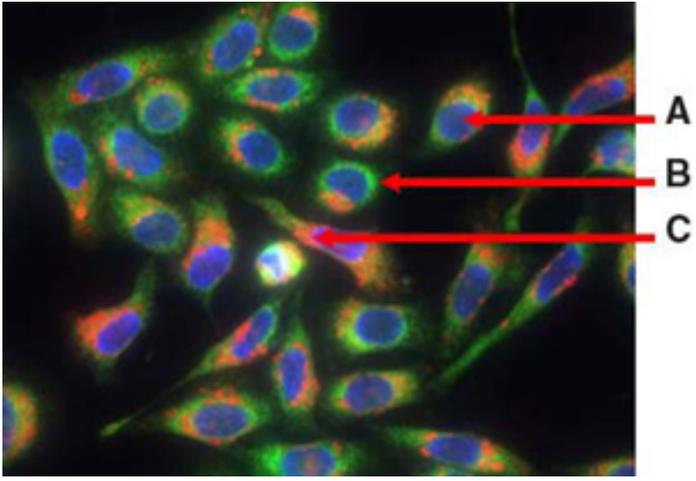


Figure 1: Live cells stained with ORGANELLE-ID-RGB[®] reagent I showing composite of (A) nuclei in blue, (B) mitochondria in green, and (C) lysosomes in red.

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

Quality Control

A sample from the ORGANELLE-ID-RGB® reagent I is used to stain HeLa cells using the procedures described in the Specific Protocol. Cells stained with ORGANELLE-ID-RGB® reagent I showing composite nuclei in blue, mitochondria in green and lysosomes in red.

Quantity

200µl (for 1000 microscopy assays).

Technical Info / Product Notes

Wavelength Maxima:

Lysosomal (Red): Excitation: 568nm; Emission: 667nm

Mitochondrial (Green): Excitation: 460nm; Emission: 560nm

Nuclear (Blue): Excitation: 350nm; Emission: 461nm

The ORGANELLE-ID-RGB® Reagent I 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 cell analysis applications involving confocal microscopy, flow cytometry, microplate readers and HCS/HTS, where consistency and reproducibility are required.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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