

ORGANELLE-ID-RGB[®] reagent II

Mixture of fluorescent dyes for detection of lysosomes, endoplasmic reticulum and nucleic acids

The ORGANELLE-ID-RGB[®] reagent II is a mixture of cell-permeable red fluorescent lysosomal dye, green fluorescent endoplasmic reticulum 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.

Ordering Information

[Order Online »](#)

ENZ-53008-C200	200µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

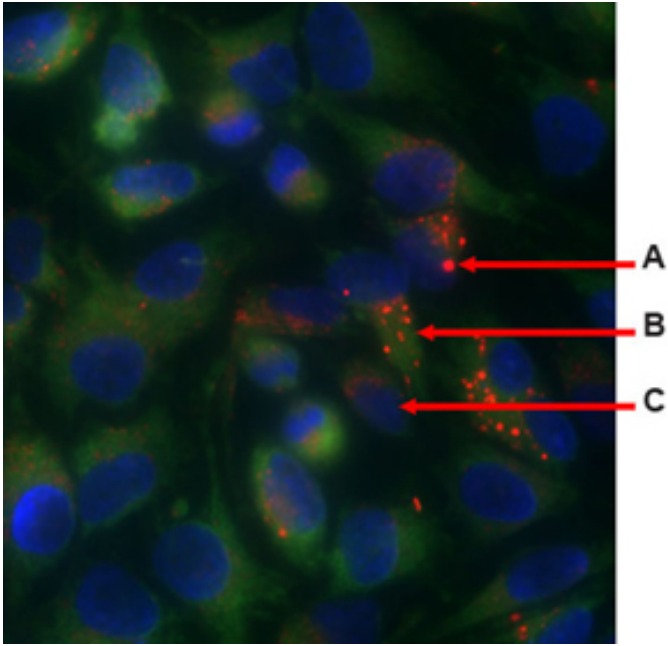


Figure 1: Live HeLa cells treated with Chloroquine and stained with ORGANELLE-ID-RGB[®] reagent II showing composite of (A) lysosomes in red, (B) endoplasmic reticula in green, and (C) nuclei in blue.

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

Purity	≥93% (HPLC)
Quantity	200µl (for 1000 microscopy assays).

Technical Info / Product Notes

Wavelength Maxima:

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

Endoplasmic Reticulum (Green): Excitation: 440nm
Emission: 565nm

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

The ORGANELLE-ID-RGB[®] Reagent II 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](mailto:info-usa@enzolifesciences.com)

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

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