

NUCLEAR-ID®

Blue/Red cell viability reagent (GFP-CERTIFIED®)

Distinguishes between live vs dead cells in a single reagent.

The NUCLEAR-ID® Blue/Red cell viability reagent (GFP-CERTIFIED®) (Prod. No. ENZ-53005) is a mixture of a blue fluorescent cell-permeable nucleic acid dye and a red fluorescent cell-impermeable nucleic acid dye that is suited for staining dead nuclei. The staining pattern arising from the simultaneous combination of these two dyes permits determination of live and dead cell populations by fluorescence/confocal microscopy. The reagent, supplied as a 1000x solution, is sufficient for 1000 microscopy assays. The single-tube format makes this cell viability reagent easy to use. It leaves the cytoplasm unstained, potentially allowing visualization of other cell markers. This reagent is specifically designed for use with GFP-expressing cell lines.

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-53005-C100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

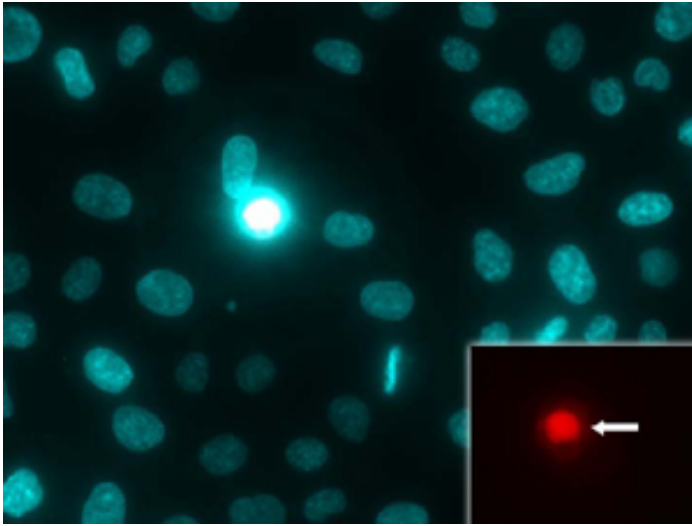


Figure 1: NUCLEAR-ID[®] Blue/Red dye is detected as blue-stained nuclei in live cells and fluorescent-red nuclei in dead cells (inset, arrow).

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	Flow Cytometry, Fluorescence microscopy, Fluorescent detection
Purity	≥93% (HPLC)
Quantity	100µl (for 1000 microscopy assays).

Technical Info / Product Notes

Application Notes:

[Use of 3D Cultured Human iPSC-Derived Hepatocytes for Long-Term Hepatotoxicity Studies](#)

[Image-Based Analysis of a Human Neurosphere Stem Cell Model for the Evaluation of Potential Neurotoxicants](#)

[Validation of a Novel Tumoroid-Based Cell Culture Model to Perform 3D in vitro Cell Signaling Analyses](#)

Wavelength Maxima

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

Dead (Red): Excitation: 571, 619nm; Emission: 639nm

NUCLEAR-ID[®] Blue/Red cell viability reagent 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.



Last modified: May 29, 2024

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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