

NUCLEAR-ID[®]

Red/Green cell viability reagent

Distinguishes between live vs dead cells in a single reagent

The NUCLEAR-ID[®] Red/Green cell viability reagent (Prod. No. ENZ-53006) is a mixture of a red fluorescent cell-permeable nucleic acid dye and a green 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.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-53006-C100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

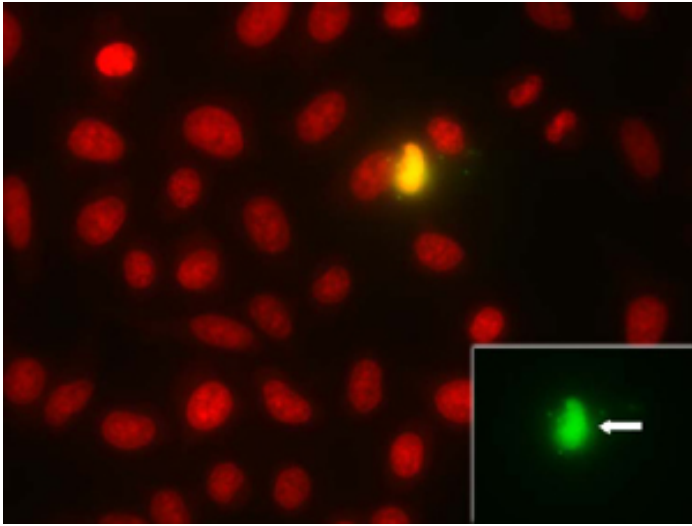


Figure 1: NUCLEAR-ID[®] Red/Green dye is detected as red-stained nuclei in live cells and fluorescent-green nuclei in dead cells (inset, arrow).

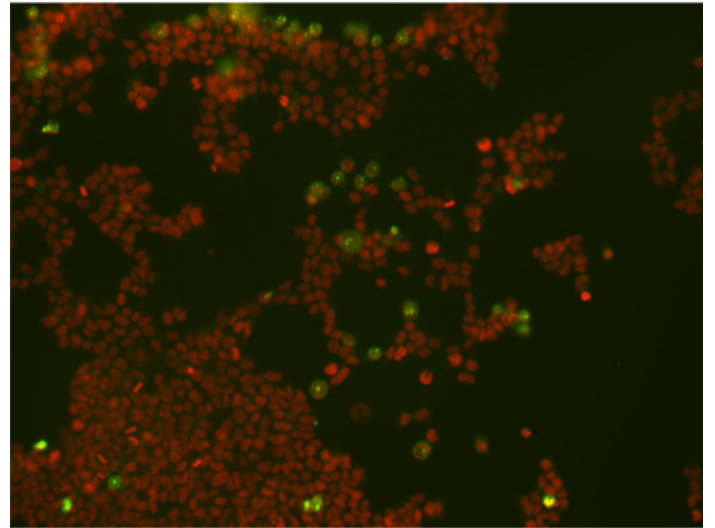


Figure 2: NUCLEAR-ID[®] Red/Green cell viability reagent (Prod. No. ENZ-53006) was used to stain monolayer cultured HeLa cell at 1:1000 dilution. The same field was imaged with appropriate filters (Red – live cell and Green – dead cell) separately and overlaid to reveal viability of cell population.

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

Wavelength Maxima

Live (Red): Excitation: 568nm; Emission: 632nm

Dead (Green): Excitation: 503nm; Emission: 524nm

The NUCLEAR-ID[®] Red/Green 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-
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)