

# NUCLEAR-ID<sup>®</sup> Blue

## DNA stain (GFP-CERTIFIED<sup>®</sup>)

### Cell permeable DNA stain that can be used for a wide range of applications

The NUCLEAR-ID<sup>®</sup> Blue DNA stain is a cell permeable dye, designed for use in a range of fluorescence detection technologies, in the discrimination of nucleated cells. It is resistant to photobleaching and is suitable for live-cell staining of nuclei. Also this dye provides a convenient approach for studying the induction and inhibition of cell cycle progression by flow cytometry. Potential applications of this reagent for live-cell studies are in the determination of cellular DNA content and cell cycle distribution, for the detection of variations in growth patterns, for monitoring apoptosis, and for evaluating tumor cell behavior and suppressor gene mechanisms.

Citations: 3

[View Online »](#)

### Ordering Information

[Order Online »](#)

ENZ-CHM103-0200	200µl
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)

- This blue fluorescent specific DNA dye is stable and has high purity
- Live, permeable and fixed cells can be analyzed
- No photobleaching effect
- No RNase treatment is required
- Multiplex with GFP and red dyes
- Validated for a wide range of cell densities
- Quick and easy to use!

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
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

Appearance	Frozen liquid.
Application	Flow Cytometry, Fluorescent detection
Emission Maximum	461nm
Excitation Maximum	350nm
Purity	≥95% (HPLC)
Quantity	200µl

Technical Info / Product Notes	The NUCLEAR-ID <sup>®</sup> Blue DNA stain is a member of the CELLESTIAL <sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL <sup>®</sup> 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.
--------------------------------	---



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-eu@enzolifesciences.com)

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