# TOTAL-NUCLEAR-ID® green/red nucleolar/nuclear detection kit

Enzo Life Sciences' TOTAL-NUCLEAR-ID® Green/Red Nucleolar/Nuclear Detection Kit contains two proprietary dyes suitable for simultaneous live-cell staining of nucleoli and nuclei. The dyes allow examination of nucleolar dynamic changes in intracellular distribution, trafficking and localization arising from biological processes such as the cell cycle and ribosome biogenesis. The kit is compatible with most fluorescence detection systems, including conventional and confocal fluorescence microscopes, as well as High Content Screening (HCS) platforms. This kit is specifically designed for visualizing nucleoli and nuclei in living cells. A control nucleolus perturbation agent, actinomycin D, is provided for monitoring changes in nucleolar dynamics. Potential applications for this kit include monitoring impaired ribosome biogenesis, inhibition transcription, cell cycle dynamics and cellular stress, as well as the distribution, trafficking and dynamics of nucleolar proteins, the distribution of viral proteins, and potentially as an aid in identifying cancer cells.

Citations: 6

View Online »

**Ordering Information** 

**Order Online** »

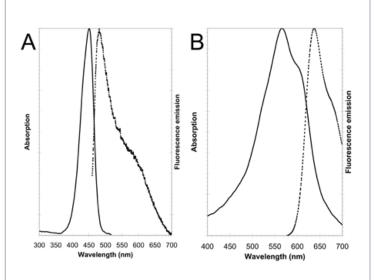
ENZ-51006-500

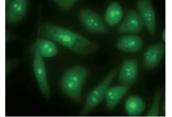
1Kit

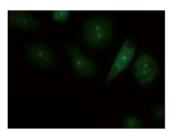
Manuals, SDS & CofA

View Online »

- Simultaneous staining of both the nucleolus and the nucleus in live cells
- High resistance to photobleaching and concentration quenching, ensuring strong, consistent fluorescence signal, even after extended viewing periods
- Validated for utility in live cell imaging applications, demonstrating appropriate response to treatment with well characterized nucleolusperturbation agents
- Stringently manufactured, to control and eliminate non-specific assay artifacts

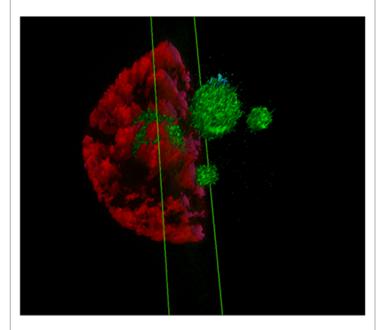


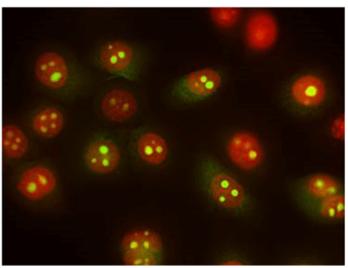




Absorption and emission spectra of A) NUCLEOLAR-ID <sup>®</sup> Green Detection Reagent and B) NUCLEAR-ID <sup>®</sup> Red Detection Reagent.

Cells stained with NUCLEOLAR-ID<sup>®</sup> Green Detection Reagent show maximal fluorescence signal within the nucleoli, and faint fluorescence throughout the nucleus and cytoplasm (A). When the cells are treated with low doses of actinomycin D, loss of nucleolar staining is observed (B). Counterstaining with NUCLEAR-ID<sup>®</sup> Red Detection Reagent facilitates highlighting the nucleoli relative to the weak cytoplasmic staining.





3D reconstruction of nucleus stained red and nucleolus stained green in U2OS cells.

Live U2OS cells, stained with TOTAL-NUCLEAR-ID<sup>®</sup> Green/Red Nucleolar/ Nuclear Detection Kit demonstrates staining of the nucleus (red) and nucleolus (yellow-green) via fluorescence microscopy. As an RNA stain, NUCLEOLAR-ID<sup>®</sup> Green dye does display some cytoplasmic staining as well, but NUCLEAR-ID<sup>®</sup> Red dye facilitates unambiguous identification of the nucleoli as green fluorescence signal within the confines of the highlighted red nuclear fluorescence signal.

# **Handling & Storage**

**Use/Stability** With proper storage, the kit components are stable up to the date noted on the product

label. Store kit at -20°C in a non-frost free freezer, or -80°C for longer term 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** Fluorescence microscopy, Fluorescent detection

Application Notes For visualizing nucleoli and nuclei simultaneously in living

cells.

Contents NUCLEOLAR-ID<sup>®</sup> Green Detection Reagent, 50 μL

NUCLEAR-ID® Red Detection Reagent, 50 µL

Actinomycin D Control, 125 μg

10X Assay Buffer, 15 mL

Quality Control A sample from each lot of TOTAL-NUCLEAR-ID<sup>®</sup> Green/

Red Nucleolus/Nucleus Detection Kit is used to stain HeLa cells using the procedures described in the user manual. Cells stained with NUCLEOLAR-ID<sup>®</sup> Green Detection Reagent show maximal fluorescence signal within the nucleoli, with only faint fluorescence throughout the nucleus and cytoplasm. NUCLEAR-ID<sup>®</sup> Red Detection Reagent maximally stains the DNA in the cell nucleus. Cells induced by actinomycin D show reduction of

nucleolar signal.

Quantity 500 assays

### **Technical Info / Product Notes**

The TOTAL-NUCLEAR-ID<sup>®</sup> green/red nucleolar/nuclear detection kit 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.

Last modified: May 29, 2024



eu@enzolifesciences.com