



Product Manual

NUCLEOLAR-ID® Green Detection Kit

Catalog #: ENZ-51009-500

500 assays



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON- INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Introduction	2
Precautions	3
Materials Supplied.....	4
Other Materials Needed	4
Reagent Preparation	5
Cell Preparations	6
Appendices	7
Results.....	8
References.....	9
Trouble Shooting Guide	10
Contact Information.....	14

INTRODUCTION

Enzo Life Sciences' NUCLEOLAR-ID® Green Detection Kit contains a proprietary dye suitable for live-cell staining of nucleoli. The dye allows examination of nucleolar dynamic changes in intracellular distribution, trafficking and localization arising from biological processes such as the cell cycle and ribosome biogenesis. Historically, nucleolus imaging approaches have required much more laborious and time consuming methods, such as fluorescently labeled RNA microinjection, fluorescence in situ hybridization (FISH) or use of fluorescent protein-tagged RNA binding proteins (GFP or YFP constructs).

The nucleolus represents a highly dynamic nuclear domain arising from an equilibrium between the level of ribosomal RNA synthesis and the efficiency of ribosomal RNA processing.^{1,2} Although the nucleolus is primarily associated with ribosome biogenesis, several lines of evidence demonstrate that it has additional functions, such as regulation of mitosis, cell-cycle progression and proliferation, and many forms of stress response and biogenesis of multiple ribonucleoprotein particles. Ribosome biogenesis is regulated throughout interphase and ceases during mitosis. Thus, there is a direct relationship between cell growth and nucleolar activities. Nucleoli are well known to be dramatically modified in cancer cells. Additionally, a large number of key proteins from both DNA- and RNA-containing viruses are localized in the nucleolus, including the HIV-1 (human immunodeficiency virus type 1) Rev and Tat proteins. Targeting of viral proteins to the nucleolus not only facilitates virus replication, but may also be required for pathogenic processes. The nucleolus is also a sensor of stress due to the redistribution of the ribosomal proteins in the nucleoplasm through its disruption.

The NUCLEOLAR-ID® Green Detection Kit is specifically designed for visualizing nucleoli in living cells by fluorescence microscopy. The dye in the kit is resistant to photobleaching, facilitating its use in imaging applications. This nucleolar stain is compatible with common live-cell nuclear stains such as Hoechst 33342, Draq5™, Vybrant® DyeCycle®Ruby and Enzo's Nuclear-ID® Red dye. The kit includes a control nucleolus perturbation agent, actinomycin D, for monitoring changes in nucleolar dynamics. Actinomycin D, an antibiotic, is a DNA-dependent RNA synthesis inhibitor. Nucleolar synthesis of ribosomal RNA is especially sensitive to actinomycin D. Potential applications for this kit include monitoring of impaired ribosome biogenesis, inhibition of transcription, cell cycle dynamics, cellular stress, distribution, trafficking and dynamics of nucleolar proteins, distribution of viral proteins, and potentially, as an aid to identify cancer cells.

SAFETY WARNINGS & PRECAUTIONS

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.



Handle
with care

- This product is for research use only and is not intended for diagnostic purposes.
- The NUCLEOLAR-ID[®] Green detection reagents contain DMSO which is readily absorbed through the skin. DMSO is harmful if ingested or absorbed through the skin and may cause irritation to the eyes. Observe appropriate precautions when handling these reagents.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas. All blood components and biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.
- To avoid photobleaching, perform all manipulations in low light environments or protected from light by other means.



Avoid
freeze /
thaw cycles

MATERIALS SUPPLIED

All reagents are shipped on dry ice. Upon receipt, the kit should be stored at $\leq -20^{\circ}\text{C}$, protected from light. When stored properly, these reagents are stable for at least twelve months. Avoid repeated freezing and thawing.

Reagents provided in the kit are sufficient for approximately 500 assays using either live, adherent cells or cells in suspension.

Reagent	Quantity
NUCLEOLAR-ID [®] Green Detection Reagent	50 μL
Actinomycin D Control	125 μg
10X Assay Buffer	15 mL



Reagents require separate storage conditions.

OTHER MATERIALS NEEDED

1. Standard fluorescence microscope
2. Calibrated, adjustable precision pipettors, preferably with disposable plastic tips
3. Adjustable speed centrifuge with swinging buckets (for suspension cultures)
4. Glass microscope slides
5. Glass cover slips
6. Deionized water
7. Anhydrous DMSO (optional)
8. Growth medium (e.g., Dulbecco's Modified Eagle Medium, D-MEM)

METHODS AND PROCEDURES

NOTE: Allow all reagents to thaw at room temperature before starting with the procedures. Upon thawing, gently hand-mix or vortex the reagents prior to use to ensure a homogenous solution. Briefly centrifuge the vials at the time of first use, as well as for all subsequent uses, to gather the contents at the bottom of the tube.

REAGENT PREPARATION

1. Positive Control

The antibiotic, actinomycin D is a DNA-dependent RNA synthesis inhibitor. Nucleolar synthesis of ribosomal RNA is especially sensitive to actinomycin D. The rearrangement of the nucleolus due to actinomycin D treatment is widely used to examine the localization of various nucleolar components, including nucleolar proteins.³ Typically, at higher doses of the drug (4-10 $\mu\text{g}/\text{mL}$ for 4 hours), the nucleolus in mammalian cells disappears, or where still present, is dramatically reduced in amount, while at lower concentrations (1-4 $\mu\text{g}/\text{mL}$ for 4 hours), less dramatic reduction in nucleolar amount is often observed.

The actinomycin D provided in the kit may be used as a positive control for reducing nucleoli size and number. It is supplied lyophilized (125 μg) and should be centrifuged briefly to gather the material at the bottom of the tube. Reconstitute the lyophilized material in 250 μL deionized water for a 0.5 mg/mL stock solution. Vortex vigorously, allow to sit for 15 minutes, then vortex again to completely bring it into solution. It is recommended that treatment with the agent be performed using 1-5 $\mu\text{g}/\text{mL}$ final concentration in order to observe changes in nucleolar morphology. Unused stock actinomycin D may be stored in small aliquots at -20°C for several weeks.

2. 1X Assay Buffer

Allow the 10X Assay Buffer to warm to room temperature. Make sure that the reagent is free of any crystallization before dilution. Prepare enough 1X Assay Buffer for the number of samples to be assayed by diluting each milliliter (mL) of the 10X Assay Buffer with 9 mL of deionized water

3. NUCLEOLAR-ID® Green Detection Reagent

The concentration of NUCLEOLAR-ID® Green dye for optimal staining will vary depending upon the application. Suggestions are provided to use as guidelines, though some modifications may be required depending upon the particular cell type employed and other factors such as the permeability of the dye to the cells or

tissues. To reduce potential artifacts from overloading of the cells, the concentration of the dye should be kept as low as possible.

Prepare sufficient amount of NUCLEOLAR-ID® Green Detection Reagent for the number of samples to be assayed as follows: To each mL of 1X Assay Buffer (see preparation in step 2, page 5) or cell culture medium containing serum, add 1 µL of NUCLEOLAR-ID® Green Detection Reagent. Serum may be included, if preferred.

NOTE: (a) If double staining is desired, the dyes may be combined into one staining solution or each may be used separately.

(b) The concentration of the NUCLEOLAR-ID® Green dye may be decreased to equal the staining intensity of the Nuclear-ID® Red dye (see Prod. No. ENZ-51006).

CELL PREPARATIONS

Cells should be maintained via standard tissue culture practices. Positive control cells should be pretreated with the actinomycin D control for 2-6 hours. Response to actinomycin D is time and concentration dependent and may also vary significantly depending upon cell type and cell line. Negative control cells should be treated with a vehicle (DMSO, media or other solvent used to reconstitute or dilute an inducer or inhibitor) for an equal length of time under similar conditions.

STAINING LIVE, ADHERENT CELLS

1. Grow cells on cover slips inside a Petri dish filled with the appropriate culture medium. When the cells have reached the desired level of confluence, carefully remove the medium.
2. Dispense sufficient volume of NUCLEOLAR-ID® Green Detection Reagent (see in step 3, page 5) to cover the monolayer cells (~100 µL of labeling solution for cells grown on an 18 X 18 mm coverslip).
3. Protect samples from light and incubate for 15-30 minutes at 37°C.
4. Wash the cells with 100 µL 1X Assay Buffer. Remove excess buffer and place coverslip on slide.
5. Analyze the stained cells by wide-field fluorescence or confocal microscopy (60X magnification recommended). Use a standard FITC filter set for imaging the nucleolus, and a standard Rhodamine or Texas Red filter set for imaging the nucleus, if Nuclear-ID® Red Detection Reagent (see Prod. No. ENZ-51006) is also used.

STAINING LIVE CELLS GROWN IN SUSPENSION

1. Centrifuge cells for 5 minutes at 400 x g at room temperature (RT) to obtain a cell pellet.
2. Carefully remove the supernatant by aspiration and dispense sufficient volume of NUCLEOLAR-ID® Green Detection Reagent (see in step 3, page 5) to cover the dispersed cell pellet.
3. Protect samples from light and incubate for 15 to 30 minutes at 37°C.
4. Wash the cells with 100 µL 1X Assay Buffer. Remove excess buffer. Resuspend cells in 100 µL 1X Assay Buffer, then apply the cells to a glass slide and overlay with a coverslip.
5. Analyze the stained cells by wide-field fluorescence or confocal microscopy (60X magnification recommended). Use a standard FITC filter set for imaging the nucleolus, and a standard Rhodamine or Texas Red filter set for imaging the nucleus, if Nuclear-ID® Red Detection Reagent (see Prod. No. ENZ-51006) is also used.

APPENDICES

FILTER SET SELECTION

The selection of optimal filter sets for a fluorescence microscopy application requires matching the optical filter specifications to the spectral characteristics of the dyes employed in the analysis. Consult the microscope or filter set manufacturer for assistance in selecting optimal filter sets for your microscope.

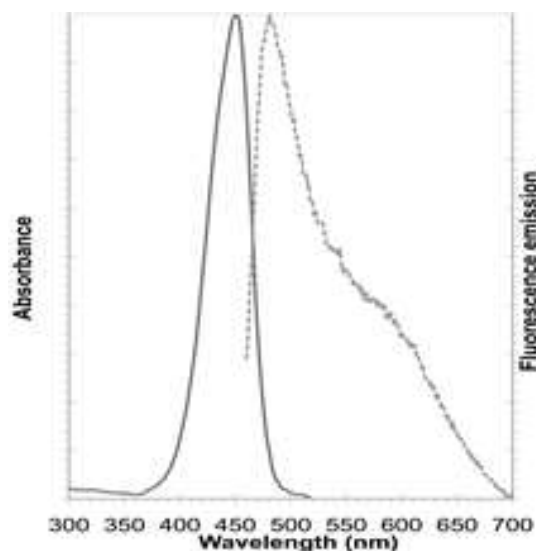


Figure 1. Absorption and fluorescence emission spectra for NUCLEOLAR-ID® Green dye. Spectrum was determined in 1X Assay Buffer.

RESULTS

Ribosomal RNAs (rRNAs) are synthesized, processed and assembled with ribosomal proteins in the nucleolus. In mammalian cells, the nucleolus is disorganized in prophase and reassembled at the end of mitosis using the nucleolar machineries from the previous cell cycle. Ribosomal DNA (rDNA) transcription is maximal in the S and G2 phases, silent in mitosis, and slowly recovers in the G1 phase of the cell cycle. The nucleolus is a prominent nuclear structure in cycling cells but of limited size in the terminal stages of cell differentiation.

Cells stained with NUCLEOLAR-ID® Green Detection Reagent show maximal fluorescence signal within the nucleoli, and faint fluorescence throughout the nucleus. Weak fluorescence is also observed throughout the cytoplasm, predominantly associated with mitochondria. The dye displays high cellular plasma and nuclear membrane permeability, and is well tolerated by living cells.

The number of nucleoli in different cell types observed with the NUCLEOLAR-ID® Green Detection Reagent will vary and they will be of different sizes as well. There appears to be an inverse relationship between size and number of nucleoli in mammalian cells.² For example, HeLa human cervical carcinoma cells, stained using the Nuclear-ID® Green Detection Reagent typically display two prominent nucleoli per cell, while U2OS human bone osteosarcoma epithelial cells are observed to contain a half dozen smaller nucleoli. Ribosomal DNA (rDNA) transcriptional arrest, induced by low doses of actinomycin D, results in loss of nucleolar staining in both cell lines as observed with the NUCLEOLAR-ID® Green Detection Reagent in the kit. The dissipation of nucleolar signal induced by actinomycin D is likely to coincide with rRNA degradation events known to occur during treatment with this drug. However, it is not definitively established whether the dye interacts with rRNA, with arginine/lysine-rich sequences in nucleolar proteins or with some other structural feature of nucleoli. NUCLEOLAR-ID® Green dye does bind nucleic acids in solution, but does not show significant selectivity towards RNA relative to DNA in such in vitro experiments.

REFERENCES

1. Hernandez-Verdun and Roussel (2003) "Regulators of nucleolar functions" *Progress in Cell Cycle Research*, Vol. 5, (Meijer, Jézéquel and Roberge, eds.) Chapter 31 pp 301-308.
2. Sirri, Urcuqui-Inchima, Roussel and Hernandez-Verdun (2008) "Nucleolus: the fascinating nuclear body" *Histochem Cell Biol* 129: 13–31.
3. Smetanaa, Buscha, Chana, Smetana and Busch (2001) "Immunocytochemical localization of nucleophosmin and RH-II/Gu. protein in nucleoli of HeLa cells after treatment with actinomycin D" *Acta Histochemica* 103, (3): 325-334

TROUBLESHOOTING GUIDE

Problem	Potential Cause	Suggestion
Nucleoli are not sufficiently stained.	Very low concentration of NUCLEOLAR-ID® Green dye was used or dye was incubated with the cells for an insufficient length of time.	Either increase the labeling concentration or increase the time allowed for the dye to accumulate in the nucleoli.
NUCLEOLAR-ID® Green dye fails to stain the nucleoli in fixed and/or permeabilized cells.	The NUCLEOLAR-ID® Green Detection Reagent provided in the kit is only suitable for live-cell staining.	Use the dye only for live-cell analysis.
Precipitate is seen in the 10X Assay Buffer.	Precipitate forms at low temperatures.	Allow solution to warm to room temperature or 37°C, then vortex to dissolve all precipitate.
The NUCLEOLAR-ID® Green stain is too bright compared to the red nuclear stain.	Different microscopes, cameras and filters may make some signals appear very bright.	Reduce the concentration of the NUCLEOLAR-ID® Green dye or shorten the exposure time.
Cells do not appear healthy.	Some cells require serum to remain healthy.	Add serum to the staining solution. Serum does not affect staining. Normal amounts of serum added range from 2% to 10%.
Actinomycin D Control does not go into solution.	The recommended volume to dissolve actinomycin D is near the limit of solubility. Ambient temperature may affect solubility.	Warm the solution to 37°C to dissolve, or dissolve in a larger volume.
Actinomycin D treated cells appear dead or are no longer attached to the surface.	The EC ₅₀ of actinomycin D may be different with different cell lines.	Lower the dose of actinomycin D, or shorten the time of exposure.



Product Manual

NOTES



Product Manual

NOTES



Product Manual

NOTES



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com

Catalog Number: ENZ-51009

Rev.: 01/23/18