

NUCLEAR-ID[®] Green

chromatin condensation detection kit

The NUCLEAR-ID[®] Green Chromatin Condensation Detection Kit detects nuclear condensation using a DNA intercalating dye that brightly stains the condensed chromatin of apoptotic cells, but only dimly stains the uncondensed chromatin of healthy living cells.

Mechanism of Action

The cell-permeable dye used in this application is an aromatic, planar cationic structure that inserts between stacked base pairs on the DNA duplex, providing an environmentally-dependent fluorescence enhancement of the dye molecules and large increases in fluorescence signal relative to the free dye in solution. Since the signal enhancement provides a proportional response, direct quantitation of DNA is possible. Further signal increase is observed upon DNA condensation during apoptosis. Considering the general mutagenic effect of nucleic acid-binding dyes, careful storage and handling of this dye is recommended.

A control apoptosis-inducing agent, staurosporine, is provided for monitoring apoptotic changes in nuclear organization. Potential applications for live-cell studies using the kit include monitoring the stages of chromatin condensation and rapid testing of compounds that induce apoptosis.

Apoptosis is recognized as a pathway of highly orchestrated signaling events, and of critical importance in biological processes and pathologies including development, aging and cancer. Nuclear condensation is one of the more prominent hallmarks of the many morphological features associated with apoptosis, including cell membrane blebbing, cell shrinkage, nucleosomal fragmentation, and formation of fragmented apoptotic bodies.

Citations: 12

[View Online »](#)

Ordering Information

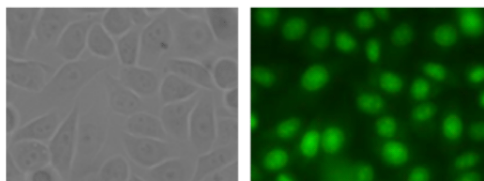
[Order Online »](#)

ENZ-51021-K200

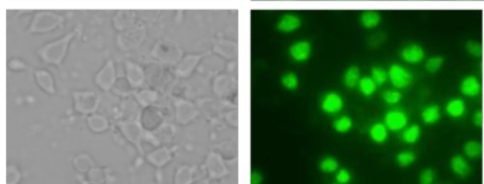
1Kit

- Dye excitable with standard 488nm laser
- Intercalating dye with superior permeability with any live cell line
- Easy no-wash mix and read protocol
- Eliminates need for specialized 350nm UV laser required for Hoechst dyes and reduces chances for channel interference
- No interference from small molecule fluorescence or cell autofluorescence

Control



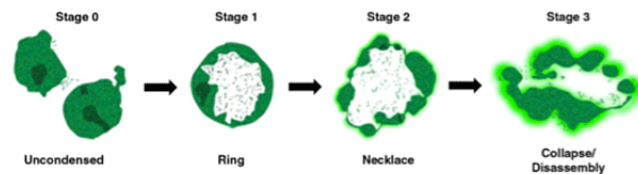
Staurosporine



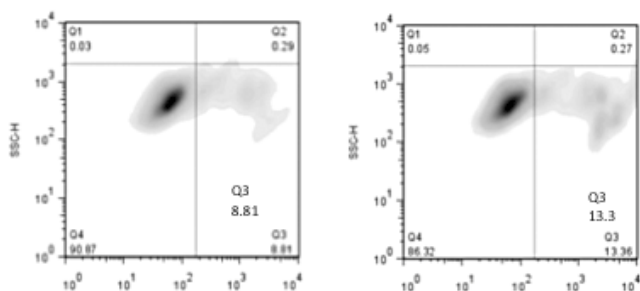
Bright Field

FITC Filter

Flow cytometry analysis of chromatin condensation in response to application of environmentally toxic compounds. Analysis was performed following 16-hour treatment of 1×10^6 cells/ml Jurkat cells with various compounds at the indicated concentration. The 488nm excitable green-emitting dye eliminates the need for specialized UV lasers required by Hoechst dye-based assays.

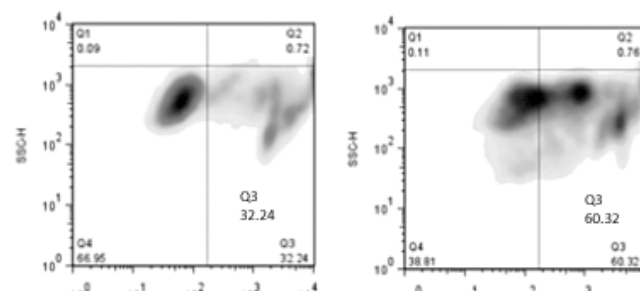


Chromatin condensation as observed by fluorescence microscopy using a standard 488nm laser. HeLa cells were treated for 4 hours with DMSO (Control) or $2 \mu\text{M}$ Staurosporine on a slide and stained with $5 \mu\text{M}$ NUCLEAR-ID[®] Green dye. The intercalating dye exhibits increased fluorescence upon chromatin condensation, a hallmark of apoptosis.



Control

10 μM BPA



10 μM Permatherin

10 μM Rotenone

Flow cytometry analysis of chromatin condensation in response to application of environmentally toxic compounds. Analysis was performed following 16-hour treatment of 1×10^6 cells/ml Jurkat cells with various compounds at the indicated concentration. The 488nm excitable green-emitting dye eliminates the need for specialized UV lasers required by Hoechst dye-based assays.

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

Application Notes This kit provides a rapid assay for nuclear condensation, a prominent hallmark of apoptosis.

Contents NUCLEAR-ID® Green Cell Cycle Detection Reagent, 100 µl Apoptosis inducer (Staurosporine), 50nmol 10X Assay Buffer, 30 ml

Quality Control Absorption peak of NUCLEAR-ID® Green dye: $\lambda_{\text{max}} = 507 \pm 4 \text{ nm}$

% purity of NUCLEAR-ID® Green dye by HPLC: $\geq 93\%$

A sample from each lot of NUCLEAR-ID® Green Cell Cycle Analysis Kit is used to analyze Jurkat cells using the procedures described in the user manual. The following results are obtained:

(a) Untreated control cells: Viable cells: $>85\%$; Apoptotic cell population: $<15\%$

(b) Apoptosis inducer treated cells: Viable cells: $>60\%$; Apoptotic cell population: $<30\%$

(c) Mean fluorescence of apoptotic cells / mean fluorescence of viable cells: >40

Quantity 200 assays



ENZO LIFE SCIENCES,
INC.
May 29, 2024
info-
usa@enzolifesciences.com

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

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