

APO-BRDU™ kit

TUNEL assay for the detection of apoptosis by flow cytometry

APO-BRDU™ Kit is a two color TUNEL (Terminal deoxynucleotide transferase dUTP Nick End Labeling) assay for labeling DNA breaks and total cellular DNA to detect apoptotic cells by flow cytometry.

The kit contains the instructions and reagents required for measuring apoptosis in cells including: positive and negative control cells for assessing reagent performance; washing, reaction and rinsing buffers for processing individual steps in the assay; terminal deoxynucleotidyl transferase enzyme (TdT), bromodeoxyuridine triphosphate (BrdUTP), and fluorescein labeled anti-BrdU antibody for labeling DNA breaks and propidium iodide/RNase A solution for counter staining the total DNA.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-0420	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application

Flow Cytometry

Contents

positive control cells, negative control cells, wash buffer, reaction buffer, TdT enzyme, Br-dUTP, rinse buffer, anti-BrdU mAb – FITC, PI/RNase staining buffer.

Quantity

50 assays.

Technical Info / Product Notes

Br-dUTP is more readily incorporated into the genome of apoptotic cells than are larger ligands like fluorescein, biotin, etc. This greater incorporation coupled with the use of a fluorescein labeled anti-BrdU monoclonal antibody (Prod. No. ALX-804-197L) gives rise to a stronger signal in flow cytometry.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
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