

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-850-042-KI01	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.



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)