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 (Br\(\text{DdUTP}\)), 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 NotesBr-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



eu@enzolifesciences.com