

### TUNEL Assay Kit - Edu-Orange ab252888

[2 Images](#)

#### Overview

|                            |  |
|----------------------------|--|
| <b>Product name</b>        | TUNEL Assay Kit - Edu-Orange   |
| <b>Detection method</b>    | Flow cytometry-fluorescent   |
| <b>Sample type</b>         | Tissue, Adherent cells, Suspension cells   |
| <b>Assay duration</b>      | Multiple steps standard assay  |
| <b>Product overview</b>    | <p>TUNEL Assay Kit - Edu-Orange ab252888 uses an EdU click chemistry method to detect fragmented DNA. The kit contains sufficient reagents to detect total/fragmented DNA in apoptotic cells in a 1 X 96-well plate or on 50 cover slips.</p> <p>The TUNEL assay is used to detect DNA fragmentation, such as in apoptosis. It uses terminal deoxynucleotidyl transferase (TdT) to catalyze the incorporation of deoxynucleotides at the free 3'-hydroxyl ends of fragmented DNA. The deoxynucleotides are then labeled in a variety of ways for detection of the degree of DNA fragmentation.</p> <p>This TUNEL assay protocol uses modified EdUTP nucleotides. A click reaction is then used to attach an orange dye (Ex 490 / Em 580, FL2 channel) to the EdUTP for detection by either flow cytometry or fluorescence microscopy. A green DNA staining dye is used for contrast (Ex 440 / Em 540)</p> <p>TUNEL assay protocol summary:</p> <ul style="list-style-type: none"> <li>- fix cells for 15 mins at room temp</li> <li>- wash cells</li> <li>- incubate with permeabilization buffer for 10 mins and wash twice</li> <li>- add TUNEL buffer and incubate for 10 mins</li> <li>- spin down cells and remove buffer</li> <li>- add TUNEL reaction cocktail and incubate for 1hr at 37°C</li> <li>- spin down cells and remove cocktail, and wash</li> <li>- add Click reaction cocktail, incubate for 30 mins and wash 3 times</li> <li>- add DNA stain, incubate for 20 mins and wash</li> <li>- analyze with flow cytometer or fluorescence microscope</li> </ul> |
| <b>Notes</b>               | <p>This product is manufactured by BioVision, an Abcam company and was previously called K191 EZClick™ TUNEL – in situ DNA Fragmentation/Apoptosis Assay Kit. K191-100 is the same size as the 100 test size of ab252888.</p>  |
| <b>Tested applications</b> | <b>Suitable for:</b> Flow Cyt, FM  |

**Platform**

Flow cytometer, Fluorescence microscope

**Properties****Storage instructions**

Please refer to protocols.

| Components                  | 100 tests |
|-----------------------------|-----------|
| 1000X Total DNA Stain       | 1 x 10µl  |
| 100X Copper Reagent         | 1 x 100µl |
| 10X Permeabilization Buffer | 1 x 25ml  |
| 10X TUNEL Reaction Buffer   | 1 x 1ml   |
| 10X Wash Buffer IV          | 1 x 25ml  |
| 20X Reducing Agent          | 1 x 500µl |
| 50X EdUTP DNA Label         | 1 x 100µl |
| Fixative Solution I         | 1 x 10ml  |
| 100X Fluorescent Azide I    | 1 x 100µl |
| TUNEL Enzyme                | 1 vial    |
| TUNEL Enzyme Buffer         | 1 x 500µl |

**Relevance**

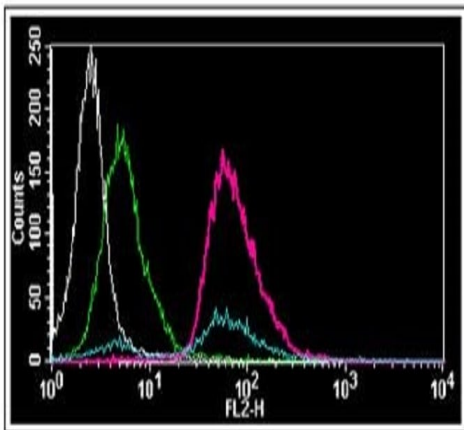
Internucleosomal DNA fragmentation is a hallmark of apoptosis in mammalian cells.

**Applications****The Abpromise guarantee**Our **Abpromise guarantee** covers the use of ab252888 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

| Application | Abreviews | Notes                                    |
|-------------|-----------|--|
| Flow Cyt    |           | Use at an assay dependent concentration. |
| FM          |           | Use at an assay dependent concentration. |

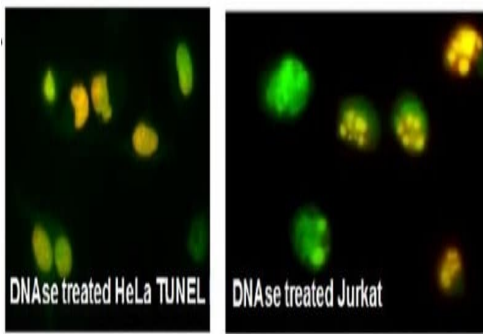
**Images**



Example Data

Detection of TUNEL-positive apoptotic strand breaks. Jurkat (Human T cell leukemia cell line from peripheral blood) cells-DNase treated ( $10^6$  cells/ml) induced strand breaks.

Unstained cells w/vehicle (white), background control cells processed for click reaction (green), negative control (untreated cells, TUNEL and click reaction; blue), DNase-treated cells (pink).



Example data

**Left panel:** DNase treated HeLa (Human epithelial cell line from cervix adenocarcinoma) cells ( $10^5$  cells/ ml). DNA staining and TUNEL and click reactions were performed.

Green: nuclear stain; Red (TUNEL positive); orange: apoptotic cells.

**Right panel:** DNase treated Jurkat (Human T cell leukemia cell line from peripheral blood) cells ( $10^6$  cells/ ml). DNA staining and TUNEL and click reactions were performed.

Green: nuclear stain; Red (TUNEL positive); orange: apoptotic cells.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors