

TUNEL Andy Fluor™ 594 Apoptosis Detection Kit

Catalog Number: A051

Table 1. Kit Components and Storage

Material	Amount	Concentration	Storage	Stability
TdT reaction buffer (Component A)	8 mL	1X solution	<ul style="list-style-type: none"> • -20 °C, • Protect from light 	<p>The product is stable for 1 year when stored as directed.</p>
TdT enzyme (Component B)	100 µL	15U/µL		
Biotin-11-dUTP (Component C)	50 µL	50X solution		
Andy Fluor 594-Streptavidin (Component D)	50 µL	100X solution		
DNase I (Component E)	10 µL	2 U/µL		
DNase I buffer (Component F)	1 mL	1X solution		
Proteinase K (Component G)	50 µL	50X solution		

Number of assays: 50 assays.

Approximate fluorescence excitation/emission maxima: Andy Fluor 594: 590/615 nm.

Introduction

Internucleosomal cleavage of DNA is a hallmark of apoptosis. DNA cleavage in apoptotic cells can be detected in situ in fixed cells or tissue sections using the terminal deoxynucleotidyl transferase (TdT) mediated dUTP nick-end labeling (TUNEL) method. TUNEL is highly selective for the detection of apoptotic cells but not necrotic cells or cells with DNA strand breaks resulting from irradiation or drug treatment.

In the TUNEL assay, TdT enzyme catalyzes the addition of labeled dUTP to the 3' ends of cleaved DNA fragments. Hapten-tagged dUTP (e.g. digoxigenin-dUTP or biotin-dUTP) can be detected using secondary reagents (e.g. anti-digoxigenin antibodies or streptavidin) for fluorescence or colorimetric detection. Alternatively, fluorescent dye-conjugated dUTP can be used for direct detection of fragmented DNA by fluorescence microscopy or flow cytometry. The TUNEL Andy Fluor™ 594 Apoptosis Detection Kit contains dUTP conjugated to biotin and Streptavidin conjugated to bright and photostable Andy Fluor™ 594 red fluorescent dye, for bright fluorescent TUNEL staining.

Materials Required but Not Provided

- Phosphate buffered saline (PBS)
- 4% formaldehyde in PBS
- 0.2% Triton X-100 in PBS
- 3% BSA in PBS
- Staining buffer: 0.6 M NaCl, 60 mM sodium citrate, 0.1% Triton X-100, 1% BSA, pH 7.4
- Hoechst 33342 counterstain (Cat No. C005)
- Antifade mounting medium
- Deparaffinization solvents (Optional)

Experimental Protocols

Sample Preparation

1. Preparation of cultured cells or fresh-frozen tissue sections

Note: Apoptotic cells can detach from adherent cell cultures and be lost during wash steps. Culture supernatants may be stained using suspension cell protocols to detect detached apoptotic cells.

- 1.1 Wash cells or sections twice in PBS.
- 1.2 Fix cells or tissues in 4% formaldehyde in PBS (pH 7.4) for 30 minutes at 4°C (not required for fixed-frozen sections)
- 1.3 Wash twice in PBS.
- 1.4 Permeabilize in PBS containing 0.2% Triton X-100 for 30 minutes at room temperature.
- 1.5 Wash twice in PBS.

2. Preparation of paraffin tissue sections

- 2.1 Deparaffinize and rehydrate tissue sections in Coplin jars at room temperature according to Tabel 2, below.

Table 2. Tissue deparaffinization procedure.

Xylenes	Xylenes	100% EtOH	100% EtOH	95% EtOH	85% EtOH	75% EtOH	1X PBS	1X PBS
5 min	5 min	5 min	5 min	5 min	3 min	3 min	5 min	5 min

- 2.2 Prepare a 1X Proteinase K solution by diluting 50X Proteinase K solution (Component G) at 1:50 in PBS. After use, aliquot any remaining stock solution and store at -20°C.
- 2.3 Permeabilize sections with 50 µL of 1X proteinase K solution for 30 minutes at room temperature. Proteinase K incubation time and temperature may require optimization depending on tissue type. Alternatively, microwave antigen retrieval protocols may be used at this step.
- 2.4 Rinse in PBS. Wash 2 x 5 minutes in PBS.

Positive Control Preparation

Note: The DNase I generates strand breaks in the DNA to provide a positive TUNEL reaction.

- 3.1 Wash sample with deionized water.
- 3.2 Incubate samples with 50 µL DNase I buffer (Component F) for 10 minutes.
- 3.3 Prepare DNase I solution according to Table 3 and mix well.

Note: Do not vortex the DNase I solution as DNase I denatures with vigorous mixing.

Table 3. DNase I solution.

Reaction Components	Number of coverslips		
	1	2	3
DNase I (Component E)	1 µL	2 µL	3 µL
DNase I buffer (Component F)	49 µL	98 µL	147 µL
Total Volume	50 µL	100 µL	150 µL

- 3.4 Remove DNase I buffer and add 50 µL of the DNase I solution to each sample and incubate for 30 minutes at room temperature.
- 3.5 Wash sample once with deionized water.

TUNEL Reaction

4.1 Incubate samples with 100 μL TdT reaction buffer (Component A) for 10 minutes.

4.2 Immediately before use, prepare the TdT reaction cocktail according to Table 4.

Table 4. TdT reaction cocktails.

Reaction Components	Number of samples				
	1	2	4	5	10
TdT reaction buffer (Component A)	47 μL	94 μL	188 μL	376 μL	470 μL
TdT enzyme (Component B)	2 μL	4 μL	8 μL	16 μL	20 μL
Biotin-11-dUTP (Component C)	1 μL	2 μL	4 μL	8 μL	10 μL
Total Volume	50 μL	100 μL	200 μL	400 μL	500 μL

4.3 Remove TdT reaction buffer and add 50 μL TdT reaction cocktail to each sample, and allow the solution to spread completely over the surface.

- a) For adherent cells or tissue sections, cover sample with a Parafilm coverslip to spread buffer evenly over cells or tissue section.
- b) For negative control samples, add 50 μL TdT reaction cocktail without TdT Enzyme.

4.4 For cell staining, incubate for 60 minutes at 37°C, protected from light. Tissue staining may require 2 hours incubation at 37°C.

- a) For adherent cells or tissue sections, perform incubation in a humid chamber.
- b) For cells in suspension, perform incubation in a microplate on a rocking platform, or resuspend cells in reaction buffer every 15 minutes by gently flicking tubes.

4.5 Wash samples 3 x 5 minutes in 3%BSA in PBS.

4.6 Prepare the Andy Fluor™ 594-Streptavidin staining solution according to Table 5.

Table 5. Andy Fluor™ 594-Streptavidin staining solution.

Reaction Components	Number of samples				
	1	2	4	5	10
Andy Fluor 594-Streptavidin (Component D)	1 μL	2 μL	4 μL	5 μL	10 μL
Staining buffer	99 μL	198 μL	396 μL	495 μL	990 μL
Total Volume	100 μL	200 μL	400 μL	500 μL	1000 μL

Staining buffer: 0.6 M NaCl, 60 mM sodium citrate, 0.1% Triton X-100, 1% BSA, pH 7.4.

4.7 Add 100 μL Andy Fluor™ 594-Streptavidin staining solution to each sample, and incubate for 30 minutes at room temperature, protected from light. Tissue staining may require 1 hours incubation at room temperature.

4.8 Wash samples 3 x 5 minutes in 3%BSA in PBS.

4.9 Counterstain samples with Hoechst 33342 if desired. Mount samples in antifade mounting medium for microscopy, or analyze cells in suspension flow cytometry.