

**VitroView™ TUNEL Apoptosis/Andy488 Kit (50 assays)**  
**Cat. No. VB-4005G**

**Introduction**

DNA fragmentation represents a characteristic hallmark of apoptosis. Terminal deoxynucleotidyl transferase dUTP nick end labeling (TUNEL) is designed to detect apoptotic cells that undergo extensive DNA degradation during the late stages of apoptosis. The method is based on the ability of terminal deoxynucleotidyl transferase (TdT) to label blunt ends of double-stranded DNA breaks independent of a template.

**Kit Components**

1. Proteinase K stock solution (20×) 0.25 mL
2. 1×Proteinase K working buffer 5 mL
3. TdT equilibration buffer 6 mL
4. TdT enzyme (20 U/μL) 20 μL
5. Biotinylated dUTP 25 μL
6. RTU Streptavidin-Andy Fluor 488 5 mL
7. TUNEL positive FFPE slides 2 slides

**Storage**

- 1×Proteinase K working buffer: Store at 4 °C.  
TdT equilibration buffer: Store at 4 °C.  
RTU Streptavidin-Andy Fluor 488: Store at 4 °C.  
TUNEL positive FFPE slides: Store at 4 °C.  
Others: Store at -20 °C.

**Protocol**

**1. Preparation of Slides**

**A. Frozen Sections**

- 1) Snap frozen fresh tissues in liquid nitrogen or isopentane pre-cooled in liquid nitrogen, embedded in OCT compound in cryomolds. Store the frozen tissue block at -80°C until ready for sectioning.
- 2) Transfer the frozen tissue block to a cryotome cryostat (e.g. -20°C) prior to sectioning and allow the temperature of the frozen tissue block to equilibrate to the temperature of the cryotome cryostat.
- 3) Section the frozen tissue block into a desired thickness (typically 5-10 μm) using the cryotome.
- 4) Place the tissue sections onto glass slides suitable for immunohistochemistry (e.g. Superfrost)
- 5) Sections can be stored in a sealed slide box at -80°C for later use.

- 6) Before staining, warm slides at room temperature for 30 minutes and fix slides by 2% neutralized formalin for 30 minutes. Rinse slides with PBS then transfer to a Coplin jar containing ice-cold 70% ethanol for 1 hour. Slides may be stored overnight in 70% ethanol at 4°C.
- 7) Wash in PBS.
- 8) Follow procedure for pretreatment as required with TdT reaction buffer.

## **B. Paraffin Sections**

- 1) Deparaffinize sections in xylene, 3×5 minutes.
- 2) Hydrate with 100% ethanol, 2×2 minutes.
- 3) Hydrate with 95% ethanol, 2×2 minutes.
- 4) Rinse in distilled water.
- 5) Prepare 1mL of 1×proteinase K working solution by mixing 50μL of 20×proteinase K solution with 950 μL of 1×Proteinase K working buffer.
- 6) Prepare carefully blot away excess water and pipette 75-100 μL of 1×proteinase K solutions to cover sections. Incubate 8-15 minutes at room temperature.
- 7) Following proteinase K treatment, wash slides 3×5 minutes with ddH<sub>2</sub>O.
- 8) (Optional) Inactivate endogenous peroxidases by covering sections with 2% hydrogen peroxide for 5 minutes at room temperature. Wash slides 3×5 minutes with ddH<sub>2</sub>O.
- 9) Follow procedure for pretreatment as required with TdT reaction buffer.

## **C. Adherent Cells**

- 1) Adherent cells may be cultured on glass chamber slides. Wash in PBS.
- 2) Fix slides by 2% neutralized formalin for 15 minutes. Rinse slides with PBS then transfer to a Coplin jar containing ice-cold 70% ethanol for 1 hour. Slides may be stored overnight in 70% ethanol at 4°C.
- 3) Wash in PBS. Inactivate endogenous peroxidases by covering sections with 2% hydrogen peroxide for 5 minutes at room temperature. Wash slides 3×5 minutes with ddH<sub>2</sub>O.
- 4) Follow procedure for pretreatment as required with TdT reaction buffer.

## **D. Suspension cells**

- 1) Harvest cells and wash them twice in PBS using centrifugation (400g for 5 minutes) to remove residual protein.
- 2) Adjust the cell concentration to 4–5×10<sup>6</sup> cells per mL in PBS.
- 3) Attach cells to slides using either the adhesion or centrifugation method.
  - a) Adhesion method
    1. Clean and label the slides.

2. Wash slides in PBS for 5 minutes at RT.
3. Place slides in a humidified box to prevent them from drying.
4. Place 20-50  $\mu\text{L}$  of the cell suspension (at least enough to cover the well) in each well of the adhesion slides and let cells adhere at room temperature (RT) for 20 minutes.
5. Fix slides by 2% neutralized formalin for 15 minutes. Rinse slides with PBS then transfer to a Coplin jar containing ice-cold 70% ethanol for 1 hour. Slides may be stored overnight in 70% ethanol at 4 °C.
6. Inactivate endogenous peroxidases by covering sections with 2% hydrogen peroxide for 5 minutes at room temperature. Wash slides 3 $\times$ 5 minutes with ddH<sub>2</sub>O.
7. Wash in PBS.
8. Follow procedure for pretreatment as required with TdT reaction buffer.

b) Centrifugation method

1. Assemble the CytoSpin centrifuge's sample chamber, filter card, slide, and racks according to the manufacturer's instructions.
2. Load 100  $\mu\text{L}$  of cells in each sample chamber.
3. Centrifuge the slides at 600 rpm for 2-4 minutes.
4. Remove the slides from the rack and place them on a staining rack.
5. Fix slides by 2% neutralized formalin for 15 minutes. Rinse slides with PBS then transfer to a Coplin jar containing ice-cold 70% ethanol for 1 hour. Slides may be stored overnight in 70% ethanol at 4 °C.
6. Wash in PBS. Inactivate endogenous peroxidases by covering sections with 2% hydrogen peroxide for 5 minutes at room temperature. Wash slides 3 $\times$ 5 minutes with ddH<sub>2</sub>O.
7. Follow procedure for pretreatment as required with TdT reaction buffer.

## 2. TUNEL Reaction

- 1) Carefully blot away excess water then cover sections with TdT equilibration buffer for 10 minutes at room temperature.
- 2) Preparation of TdT reaction buffer:

	2 samples	4 samples	10 samples
TdT equilibration buffer	24 $\mu\text{L}$	48 $\mu\text{L}$	94 $\mu\text{L}$
TdT enzyme	0.5 $\mu\text{L}$	1 $\mu\text{L}$	2.5 $\mu\text{L}$
Biotinylated dUTP	0.8 $\mu\text{L}$	1.6 $\mu\text{L}$	4 $\mu\text{L}$

Mix well. Prepare fresh from stock solution prior to use.

- 3) Remove TdT equilibration buffer and cover sections with 10  $\mu\text{L}$  of TdT reaction buffer. Incubate slides in a humidified chamber for 30 minutes at 37 °C. In order to conserve reagents a reduced volume of TdT buffer may be carefully covered with a glass coverslip during the incubation. Take care to avoid trapping air bubbles which may lead to staining artifacts.
- 4) Stop reaction by incubating slides 2 $\times$ 10 minutes in 1 $\times$ SSC.

## 3. Visualized with Andy Fluor 488

- 1) Rinse slides in PBS then block nonspecific binding by covering tissue sections with 2% BSA solution for 10 minutes at room temperature.
- 2) Rinse in PBS for 3 $\times$ 2 minutes.

- 3) **Detection:** Incubate sections with 3-4 drops of RTU streptavidin-Andy Fluor 488 for 30 minutes at room temperature.
- 4) Rinse in PBS for 3×2 minutes.
- 5) **Counterstain:** Dilute the DAPI stock solution to 300 nM in PBS. Add approximately 300 µL of this dilute DAPI staining solution to the section/cover slip, making certain that the tissue/cells are completely covered.
- 6) Incubate for 1 – 5 minutes.
- 7) Rinse the sample several times in PBS. Drain excess buffer from the section/cover slip and mount. We recommend using a mounting medium with an antifade reagent.
- 8) View the sample using a fluorescence microscope with appropriate filters.

**Negative Control:** As a negative control, omit the TdT enzyme from the TdT reaction buffer.

**Note:** This product is intended for research purposes only. This product is **not** intended to be used for therapeutic or diagnostic purposes in humans or animals.

**Precautions:** Handle with care. Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.