

## Cell Meter™ Annexin V Binding Apoptosis Assay Kit

*\*Orange Fluorescence Optimized for Flow Cytometry \**

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 22825 (100 assays)	Keep at 4 °C and avoid exposure to light	Flow cytometry Fluorescence microscope

### Introduction

Annexins are a family of calcium-dependent phospholipid-binding proteins. They are abundant in eukaryotic organisms belonging to a family of ubiquitous cytoplasmic proteins involved in signal transduction. Annexin V's preferential binding partner is phosphatidylserine (PS), which is usually kept on the inner-leaflet (the cytosolic side) of cell membranes. In apoptosis, PS is transferred to the outer leaflet of the plasma membrane. The appearance of phosphatidylserine on the cell surface is a universal indicator of the initial/intermediate stages of cell apoptosis and can be detected before morphological changes can be observed. Our Cell Meter™ assay kits are a set of tools for monitoring cell apoptosis through measuring the translocation of phosphatidylserine (PS).

This kit uses our proprietary orange fluorescent Annexin V-iFluor™ 555 PS sensor that specifically binds PS with orange fluorescence. The spectral property is almost identical to those of Cy3® or Alexa Fluor® 555, making it convenient to be used for common fluorescence instruments equipped with the light sources and filters for Cy3® or Alexa Fluor® 555 (Cy3® or Alexa Fluor® 555 are the trademarks of GE Healthcare and Invitrogen respectively). This kit provides all the essential components with an optimized protocol for flow cytometric applications.

### Kit Key Features

<b>Non-Radioactive:</b>	No special requirements for waste treatment.
<b>Convenient:</b>	All essential assay components are included.
<b>Optimized Performance:</b>	Provide optimal conditions for detecting the translocation of phosphatidylserine.
<b>Enhanced Value:</b>	Less expensive than the sum of individual components

### Kit Components

Components	Amount
Component A: Annexin V-iFluor™ 555 (100X stock solution)	1 vial (200 µL/vial)
Component B: Assay Buffer	50 mL

### Assay Protocol

#### Brief Summary

**Prepare cells with test compounds (200 µL/sample) → Add Annexin V-iFluor™ 555 assay solution  
→ Incubate at room temperature for 30-60 minutes → Analyze cells using a flow cytometer  
or Fluorescence microscope with Ex/Em = 555/ 590 nm filter set**

#### 1. Prepare and incubate cells with Annexin V-iFluor™ 555:

- 1.1 Treat cells with test compounds for a desired period of time (4-6 hours for Jurkat cells treated with staurosporine) to induce apoptosis.
- 1.2 Centrifuge the cells to get  $1-5 \times 10^5$  cells/tube.
- 1.3 Resuspend cells in 200 µL of Assay Buffer (Component B).
- 1.4 Add 2 µL of Annexin V-iFluor™ 555 (Component A) into the cells.
- 1.5 Incubate at room temperature for 30 to 60 minutes, protected from light.
- 1.6 Add 300 µL of Assay Buffer (Component B) to increase volume before analyzing the cells with a flow cytometer or fluorescence microscope (see Step 1.7 below).
- 1.7 Monitor the fluorescence intensity at Ex/Em = 555/590 nm by using a flow cytometer or a fluorescence microscope (See Step 2 or 3 below).

**2. Analyze by using a flow cytometer:**

Quantify Annexin V- iFluor™ 555 binding by using a flow cytometer at Ex/Em = 555/590 nm.

*Note: Annexin V binding flow cytometric analysis on adherent cells is not routinely tested since specific membrane damage may occur during cell detachment or harvesting. However, methods for utilizing Annexin V for flow cytometry on adherent cell types have been previously reported by Caciola-Rosen et al. and van Engeland et al (see Refs 1 and 2).*

**3. Analyze by using a fluorescence microscope:**

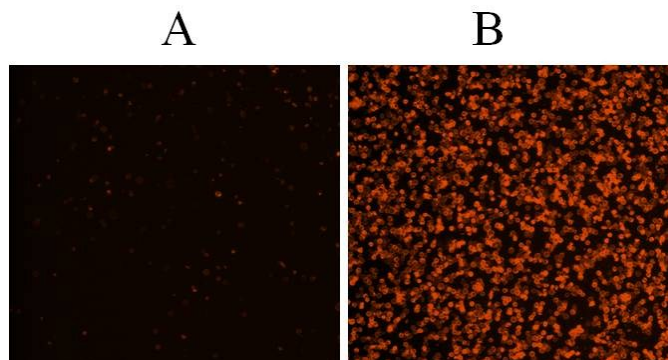
- 3.1 Pipette the cell suspension from Step 1.5, rinse 1-2 times with assay buffer, and then resuspend the cells with assay buffer. Add the cells on a glass slide that is covered with a glass cover slip.

*Note: For adherent cells, it is recommended to grow the cells directly on a cover slip. After incubation with Annexin V-iFluor™ 555 (Step 1.5), rinse 1-2 times with assay buffer, and add assay buffer back to the cover slip. Invert cover slip on a glass slide and visualize the cells. The cells can also be fixed in 2% formaldehyde after the incubation with Annexin V-iFluor™ 555 and visualized under a microscope.*

- 3.2 Analyze the apoptotic cells with Annexin V-iFluor™ 555 under a fluorescence microscope using the Violet channel. Measure the cell viability by using the Cy5 channel when Nuclear Red™ DCS1 (AAT catalog# 17552) is added into the cells. The orange staining on the plasma membrane indicates the Annexin V-iFluor™ 555 binding to PS on cell surface.

**Data Analysis**

In live non-apoptotic cells, Annexin V-iFluor™ 555 detects innate apoptosis in non-induced cells, which is typically 2-6% of all cells. In apoptotic cells, Annexin V-iFluor™ 555 binds to phosphatidylserine, which is located on the outer leaflet of the cell membrane, resulted in increased staining intensity.



**Figure1.** Images of Jurkat cells in a Costar black wall/clear bottom 96-well plate stained with the Cell Meter™ Annexin-V Binding Apoptosis Assay Kit \*Orange Fluorescence\* A: Untreated control cells B: Cells treated with 1 μM staurosporine for 5 hours.

**References**

1. van Engeland M, Ramaekers FCS, Schutte B, Reutelingsperger CPM: A novel assay to measure loss of plasma membrane asymmetry during apoptosis of adherent cells in culture. *Cytometry* 24:131–139, 1996.
2. L Caciola-Rosen, A Rosen, M Petri, and M Schlissel. Surface blebs on apoptotic cells are sites of enhanced procoagulant activity: implications for coagulation events and antigenic spread in systemic lupus erythematosus. *Proc Natl Acad Sci U S A.* 1996 February 20; 93(4): 1624–1629.
3. Hanshaw RG, Lakshmi C, Lambert TN, Johnson JR, Smith BD. (2005) Fluorescent detection of apoptotic cells by using zinc coordination complexes with a selective affinity for membrane surfaces enriched with phosphatidylserine. *Chembiochem*, 6, 2214.
4. Koulov AV, Stucker KA, Lakshmi C, Robinson JP, Smith BD. (2003) Detection of apoptotic cells using a synthetic fluorescent sensor for membrane surfaces that contain phosphatidylserine. *Cell Death Differ*, 10, 1357.

**Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest®. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.**