

## Cell Meter™ FITC-Annexin V Binding Apoptosis Assay Kit

*\*Optimized for Flow Cytometry\**

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 22839 (100 assays)	Keep at 4 °C and avoid exposure to light	Flow Cytometer

### 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 Annexin V- FITC sensor that specifically binds to PS for the early stages of apoptosis, and in pair with uptake of propidium iodide (PI) for late-stage apoptosis. This kit provides all the essential components with an optimized protocol for flow cytometric and fluorescence microscopy applications.

### Kit Key Features

<b>Non-Radioactive:</b>	No special requirements for waste treatment.
<b>Convenient:</b>	Include all essential assay components.
<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- FITC (100X stock solution)	1 vial (200 µL/vial)
Component B: Assay Buffer	50 mL
Component C: 100X Propidium Iodide	1 vial (100 µL)

### Assay Protocol

#### Brief Summary

**Prepare cells with test compounds (200 µL/sample) → Add Annexin V-FITC assay solution  
→ Incubate at room temperature for 30-60 minutes → Analyze cells with a  
flow cytometer using FL1 channel (Ex/Em = 490/525 nm)**

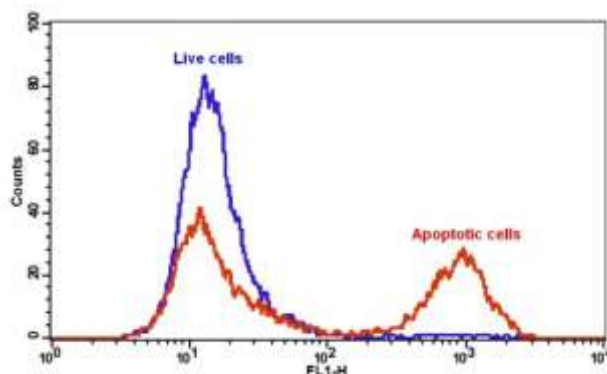
1. Treat cells with test compounds for a desired period of time (4-6 hours for Jurkat cells treated with camptothecin) to induce apoptosis.  
*Note: Annexin V flow cytometric analysis on adherent cells is not routinely tested since specific membrane damage may occur during cell detachment or harvesting. 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).*
2. Centrifuge the cells to get  $2-5 \times 10^5$  cells/tube.
3. Resuspend cells in 200 µL of Assay Buffer (Component B).
4. Add 2 µL of Annexin V-FITC (Component A) into the cells.

*Optional: Add 2  $\mu$ L of 100X Propidium Iodide (Component C) for necrosis cells.*

5. Incubate at room temperature for 30 to 60 minutes, protected from light.
6. *Optional:* add 200 to 300  $\mu$ L of Assay Buffer (Component B) to increase volume before analyzing the cells with a flow cytometer (See Step 7).
7. Monitor the fluorescence intensity of Annexin V-FITC using the FL1 channel (Ex/Em = 490/525 nm), and measure the cell viability with propidium iodide using the FL2 channel.

## Data Analysis

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



**Figure1.** The detection of binding activity of Annexin V-FITC to phosphatidylserine in Jurkat cells. Jurkat cells were treated without (Blue) or with 20  $\mu$ M camptothecin (Red) in a 37 °C, 5% CO<sub>2</sub> incubator for 4-5 hours, and then dye loaded with Annexin V-FITC for 30 minutes. The fluorescence intensity of Annexin V-FITC was measured with a FACSCalibur (Becton Dickinson, San Jose, CA) flow cytometer using the FL1 channel.

## References

1. van Engeland M, Ramaekers FCS, Schutte B, Reutelingsperger CPM: A novel assay to measure loss of plasma membrane asymmetry during apoptosis of adherant cells in culture. *Cytometry* 24:131–139, 1996.
2. L Casciola-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.
5. Hall MP, Burson KK, Huestis WH. (1998) Interactions of a vesicular stomatitis virus G protein fragment with phosphatidylserine: NMR and fluorescence studies. *BiochimBiophys Acta*, 1415, 101.
6. Saurel O, Cezanne L, Milon A, Tocanne JF, Demange P. (1998) Influence of annexin V on the structure and dynamics of phosphatidylcholine/phosphatidylserine bilayers: a fluorescence and NMR study. *Biochemistry*, 37, 1403.
7. Hanada K, Pagano RE. (1995) A Chinese hamster ovary cell mutant defective in the nonendocytic uptake of fluorescent analogs of phosphatidylserine: isolation using a cytosol acidification protocol. *J Cell Biol*, 128, 793.

**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.