

## Cell Meter™ Annexin V Binding Apoptosis Assay Kit \*Green Fluorescence Optimized for Flow Cytometry\*

Catalog number: 22824  
Unit size: 100 Tests

Component	Storage	Amount
Component A: Annexin V-iFluor™ 488 (100X stock solution)	Refrigerate (2-4 °C), Avoid Light	1 vial (200 µL/vial)
Component B: Assay Buffer (4 °C)	Refrigerate (2-4 °C)	50 mL
Component C: 100X Propidium Iodide	Freeze (<-15 °C), Avoid Light	1 vial (100 µL)

### OVERVIEW

Our Cell Meter™ assay kits are a set of tools for monitoring cell viability. There are a variety of parameters that can be used for monitoring cell viability. This particular kit is designed to monitor cell apoptosis through measuring the translocation of phosphatidylserine (PS). 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. This kit uses a fluorescent Annexin V that specifically binds PS. Annexin V conjugates have been demonstrated to selectively bind PS. This particular assay kit is optimized to monitor cell apoptosis using a flow cytometer with the FITC channel (green fluorescence).

### AT A GLANCE

#### Protocol summary

1. Prepare cells with test compounds (200 µL/sample)
2. Add Annexin V-iFluor™ 488 stock solution
3. Incubate at room temperature for 30 - 60 mintues
4. Analyze cells using flow cytometer with FL1 channel (Ex/Em = 490/525 nm)

**Important** Thaw 100X Propidium Iodide (Component C) at room temperature before starting the experiment.

### KEY PARAMETERS

Instrument:	Flow cytometer
Excitation:	490 nm
Emission:	525 nm
Instrument specification(s):	FL1 channel

### PREPARATION OF CELL SAMPLES

For guidelines on cell sample preparation, please visit  
<https://www.aatbio.com/resources/guides/cell-sample-preparation.html>

### SAMPLE EXPERIMENTAL PROTOCOL

1. Treat cells with test compounds for a desired period of time (4 - 6 hours for Jurkat cells treated with camptothecin) to induce apoptosis.

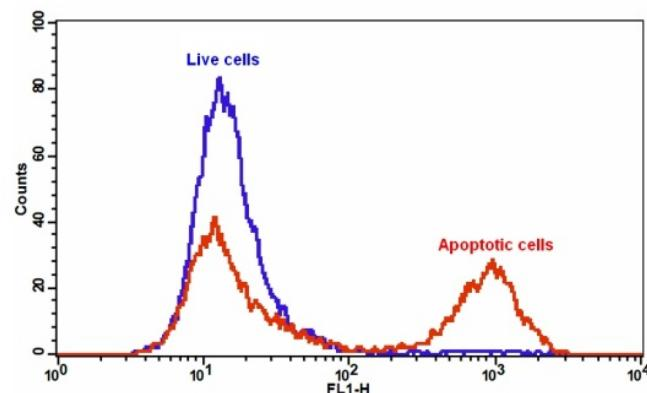
**Note** Annexin V flow cytometric analysis on adherent cells is not routinely tested since specific membrane damage may occur during cell detachment or harvesting. However, methods for utilizing Annexin V for flow cytometry on adherent cell types have been previously reported by Casiola-Rosen et al. and van Engelend et al.

2. Centrifuge the cells to get  $1 - 5 \times 10^5$  cells/tube.
3. Resuspend cells in 200 µL of Assay Buffer (Component B).

4. Add 2 µL of Annexin V-iFluor™ 488 (Component A) into the cells.
5. **Optional:** Add 2 µL of 100X Propidium Iodide (Component C) for necrosis cells.
6. Incubate at room temperature for 30 to 60 minutes, protected from light.
7. **Optional:** Add 200 to 300 µL of Assay Buffer (Component B) to increase volume before analyzing the cells with a flow cytometer.
8. Monitor the fluorescence intensity of Annexin V-iFluor™ 488 using a flow cytometer with FL1 channel (Ex/Em = 490/525 nm). Measure the cell viability using the FL2 channel when propidium iodide is added into the cells.

### EXAMPLE DATA ANALYSIS AND FIGURES

In live non-apoptotic cells, Annexin V-iFluor™ 488 detects innate apoptosis in non-induced cells, which is typically 2-6% of all cells. In apoptotic cells, Annexin V-iFluor™ 488 binds to phosphatidylserine, which is located on the outer leaflet of the cell membrane, therefore resulted in increased staining intensity



**Figure 1.** The detection of binding activity of Annexin V-iFluor™ 488 and phosphatidylserine in Jurkat cells. Jurkat cells were treated without (Blue) or with 20 µM camptothecin (Red) in a 37 °C, 5% CO<sub>2</sub> incubator for 4-5 hours, and then dye loaded with Annexin V-iFluor™ 488 for 30 minutes. The fluorescence intensity of Annexin V-iFluor™ 488 was measured with a FACSCalibur (Becton Dickinson) flow cytometer using the FL1 channel.

### DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.