

Cell Meter™ APC-Annexin V Binding Apoptosis Assay Kit *Optimized for Flow Cytometry*

Catalog number: 22837
Unit size: 100 Tests

| Component | Storage | Amount |
|--------------------------------------|-----------------------------------|-----------------|
| Component A: APC-Annexin V conjugate | Refrigerate (2-4 °C), Avoid Light | 1 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

Annexin V may be conjugated to fluorochromes including APC. This format retains its high affinity for phosphatidylserine (PS) and thus serves as a sensitive probe for flow cytometric analysis of cells that are undergoing apoptosis. Since externalization of PS occurs in the earlier stages of apoptosis, APC Annexin V staining can identify apoptosis at an earlier stage than assays based on nuclear changes such as DNA fragmentation. APC Annexin V staining precedes the loss of membrane integrity which accompanies the latest stages of cell death resulting from either apoptotic or necrotic processes. Therefore, staining with APC Annexin V is typically used in conjunction with a vital dye such as propidium iodide (PI) or 7-Amino-Actinomycin (7-AAD) to allow the investigator to identify early apoptotic cells (7-AAD negative, APC Annexin V positive). Viable cells with intact membranes exclude 7-AAD, whereas the membranes of dead and damaged cells are permeable to 7-AAD. For example, cells that are considered viable are both APC Annexin V and 7-AAD negative while cells that are in early apoptosis are APC Annexin V positive and 7-AAD negative, while cells that are in late apoptosis or already dead are both APC Annexin V and 7-AAD positive. This assay does not distinguish between cells that have undergone apoptotic death versus those that have died as a result of a necrotic pathway because in either case, the dead cells will stain with both APC Annexin V and 7-AAD. However, when apoptosis is measured over time, cells can be often tracked from APC Annexin V and 7-AAD negative (viable, or no measurable apoptosis), to APC Annexin V positive and 7-AAD negative (early apoptosis, membrane integrity is present) and finally to APC Annexin V and 7-AAD positive (end stage apoptosis and death). The movement of cells through these three stages suggests apoptosis. In contrast, a single observation indicating that cells are both APC Annexin V and 7-AAD positive, in of itself, reveals less information about the process by which the cells underwent their demise.

AT A GLANCE

Protocol summary

1. Prepare cells with test compounds (200 µL/sample)
2. Add APC-Annexin V assay solution
3. Incubate at room temperature for 20 - 60 minutes
4. Analyze cells using flow cytometer with FL4 channel (Ex/Em = 635/660 nm)

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

KEY PARAMETERS

Instrument: Flow cytometer
Excitation: FL4 channel
Emission: FL4 channel

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles.

1. APC-Annexin V stock solution (100X):
Add 200 µL PBS with 0.2% BSA into the vial of APC-Annexin V conjugate

(Component A) and mix well to make 100X APC-Annexin V stock solution.

Note Store the reconstituted 100X APC-Annexin V stock solution at 4 °C. Do Not Freeze.

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 staurosporine) 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 Engelen 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 100X APC-Annexin V stock solution into the cells.
5. **Optional:** Add 2 µL of 100X Propidium Iodide (Component C) into the cells for necrosis cells.
6. Incubate at room temperature for 20 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 APC-Annexin V using a flow cytometer with FL4 channel (Ex/Em = 635/660 nm). Measure the cell viability using FL2 channel when propidium iodide is added into the cells.

EXAMPLE DATA ANALYSIS AND FIGURES

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

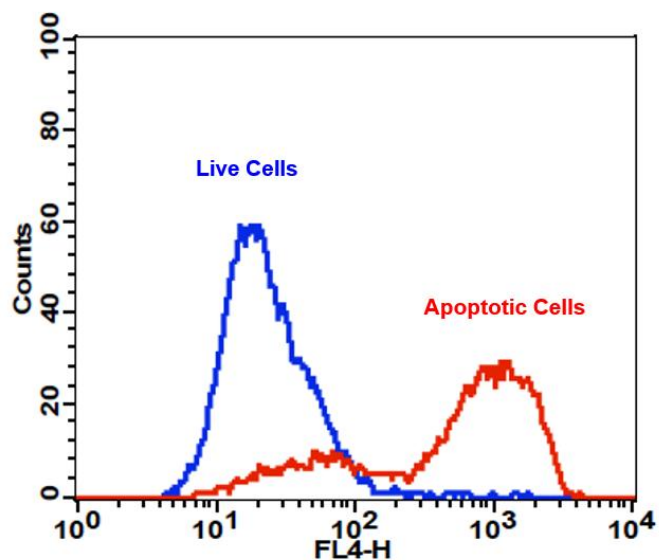


Figure 1. The detection of binding activity of APC-Annexin V to phosphatidylserine in Jurkat cells with Cell Meter™ APC-Annexin V Binding Apoptosis Assay Kit. Jurkat cells were treated without (Blue) or with 1 μ M staurosporine (Red) in a 37 °C, 5% CO₂ incubator for ~4 hours, and then dye loaded with APC-Annexin V for 30 minutes. The fluorescence intensity of APC-Annexin V was measured with a FACSCalibur (Becton Dickinson) flow cytometer using the FL4 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.