

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

Catalog number: 22837
Unit size: 100 Tests

Component	Storage	Amount (Cat No. 22837)
Component A: APC-Annexin V conjugate	Minimize light exposure, Refrigerated (2-8 °C)	1 Vial
Component B: Assay Buffer (4 °C)	Refrigerated (2-8 °C)	1 Bottle (50 mL)
Component C: 100X Propidium Iodide	Freeze (< -15 °C), Minimize light exposure	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.

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 660/20 nm filter (APC channel)

Important Note

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

KEY PARAMETERS

Flow cytometer

Emission	660/20 nm filter
Excitation	640 nm laser
Instrument specification(s)	APC channel

CELL PREPARATION

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

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

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.

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 Engeland 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 660/20 nm filter (APC channel). Measure the cell viability using 610/20 nm filter (PE-Texas Red 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.

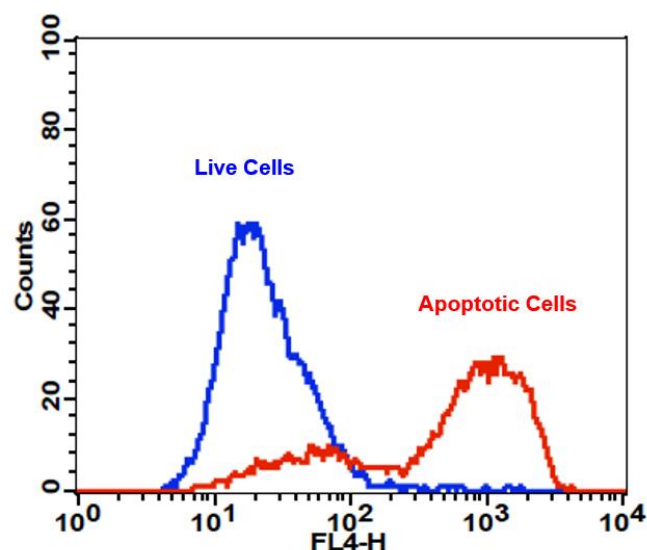


Figure 1. The detection of binding activity of APC-Annexin V to phosphatidylserine in Jurkat cells with Cell Meter™ APC-Annexin V Binding Apoptosis Assay Kit. Jurkat cells were treated without (Blue) or with 1 μ M staurosporine (Red) in a 37 °C, 5% CO₂ incubator for ~4 hours, and then dye loaded with 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.