

December 2012

Annexin V with Cell Surface Antibody Staining for Suspension Cells

Materials and Reagents

Full Name	Short Name	Catalog Number
Microwell plates (round-bottom wells) or tubes (12 x 75-mm polypropylene round-bottom test tube)		N/A
1X PBS Buffer	PBS	554781
BD Pharmingen™ stain buffer, or equivalent, 1X PBS, 2% FBS (or BSA) 0.1% NaN ₃ (pH 7.1–7.4)	Stain Buffer (FBS)	554656
Annexin V conjugate	Stain Buffer (BSA)	554657
Propidium Iodide staining solution	Annexin V	*
BD Via-Probe™ Cell Viability Solution	PI	556463
Annexin V Binding Buffer, 10X Concentrate	7-AAD	555816
N/A, not applicable	Annexin V Binding Buffer	556454
*Select at www.bdbiosciences.com		

Procedural Notes

- Investigators might want to check whether the antibody staining procedure affects cell surface levels of Annexin V by including control tubes with or without cell surface antibodies.
- Use PI with FITC-, APC-, or BD Horizon™ V450-conjugated Annexin V; use 7-AAD with PE-, Cy™ 5-, Cy™ 5.5- or BD Horizon™ V500-conjugated Annexin V.
- The following controls are used to set up compensation and quadrants:
 - Unstained cells
 - Cells stained with Annexin V-conjugate alone (no viability dye)
 - Cells stained with viability dye alone (no Annexin V-conjugate)

Procedure

- Pellet and wash cells using Stain Buffer.
- Optional: cells may be stained with antibodies against cell surface antigens prior to the Annexin V staining procedure. If you are not staining cells with antibodies, proceed to step 3.
 - Wash cells with Stain Buffer and resuspend in 100 µL in 12 x 75-mm tubes or a 96-well plate.
 - Add antibodies against cell surface antigens and incubate for 20-45 minutes in the dark (either on ice or at RT).
 - Wash cells twice (1-2 mL for tubes or 100-200 µL for 96-well plates, centrifuging at 300g) with Stain Buffer, then proceed with step 3 below.
- Wash cells twice with cold PBS and then resuspend cells in 1X Binding Buffer at a concentration of $\sim 1 \times 10^6$ cells/mL.
- Transfer 100 µL of the solution ($\sim 1 \times 10^5$ cells) to a 5-mL culture tube.
- Add Annexin V and PI or 7-AAD as described in the Technical Data Sheet or the Annexin V apoptosis detection kit manual.
- Gently mix the cells and incubate for 15 minutes at RT in the dark.



December 2012

Annexin V with Cell Surface Antibody Staining for Suspension Cells

7. Add 400 μ L of 1X Binding Buffer to each tube. Analyze by flow cytometry as soon as possible (within 1 hr).

Cy™ is a trademark of Amersham Biosciences Corp. Cy™ dyes are subject to proprietary rights of Amersham Biosciences Corp and Carnegie Mellon University and made and sold under license from Amersham Biosciences Corp. This product is licensed for sale only for research and in vitro diagnostic use. Any other use requires a commercial sublicense from Amersham Biosciences Corp, 800 Centennial Avenue, Piscataway, NJ 08855-1327, USA.



December 2012

Annexin V with Cell Surface Antibody Staining for Suspension Cells

