RayBio[®] Annexin V Apoptosis Detection Kit

(RayBright® Violet 450)

User Manual Version 1.0 Oct 31st, 2018

RayBio® Annexin V Apoptosis Detection Kit (RayBright® Violet 450)

Catalog numbers: 137-08010-25 (25 tests) 137-08010-100 (100 tests)



We Provide You With Excellent Support And Service

Tel:(Toll Free)1-888-494-8555 or 770-729-2992; Fax:770-206-2393; Web: www.raybiotech.com Email: info@raybiotech.com



RayBio® Intracellular Staining Kit

TABLE OF CONTENTS

١.	Introduction	1
II.	Kit Contents	2
III.	Assay Protocol	2
	Storage and Stability	

I. INTRODUCTION

The RayBiotech Annexin V Detection Kit is a sensitive assay for detection of cell apoptosis by flow cytometry. Annexin V is a cellular protein that specifically binds to phosphatidylserine (PS) sites on cell membrane. In live healthy cells, PS is predominantly localized along the cytosolic side of the plasma membrane. However, during induction of apoptosis or necrosis, PS translocates to the extracellular membrane. If present on the extracellular surface, the PS can be detected by fluorochrome-labeled Annexin V. With the inclusion of RayBiotech's fixable live/dead marker stain, live and dead cells can be discriminated as well. However, unlike other apoptosis assay kits using propidium iodide (PI) or 7-AAD, RayBio Annexin V apoptosis Detection Kit can be combined with immunophenotyping that requires cellular fixing of the cell populations. Combined, the RayBiotech Annexin V Detection kit allows for a fixable live/dead and apoptotic/necrotic staining kit.

II. KIT CONTENTS

Components	137-08010-25	137-08010-100	Part
	25 tests	100 tests	Number
RayBright® R780 Live (in DMSO, 1:100	25 μL	100 μL	Item A
dilution for use)			
RayBright® V450 Annexin V (5 μL/test)	125 μL	500 μL	Item B
Annexin V Binding Buffer, 5x (dilute	5 mL	20 mL	Item C
with ddH2O to 1x solution)			
Annexin V Co-staining Buffer, 5x	5 mL	20 mL	Item D
(dilute with ddH2O to 1x solution)			

- Kit can be stored at 4°C, protected from light for up to 3 months
- Make enough Live/Dead dye (Item A), 1x Binding Buffer (Item C), and Co-staining Buffer (Item D) for your assay. Return remaining stock at 4°C.

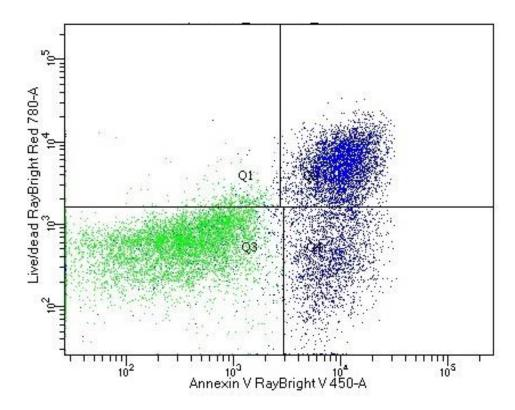
III. Assay Protocol

- 1. Transfer cells to a 96-well plate.
- 2. Wash cells twice with cold PBS, spin down cells at 300g for 10 minutes at 4°C (same centrifuge speed applies to the following steps).
- 3. Resuspend cells in 100 μ L of PBS containing diluted RayBright® 780 live/dead marker (Item A, 1:100 dilution in PBS)
- 4. Incubate for 20 minutes on ice.
- 5. Spin down the cells, wash cells twice with 200 μ L 1x Annexin V Binding Buffer (Item C)
- 6. Resuspend cells in 100 μ L of 1x annexin V binding buffer, add 5 μ L of fluorochrome-conjugated-annexin V (Item B), incubate at room temperature for 15 minutes.
- 7. Wash the cells once with 200 μ L 1x Annexin V Binding Buffer, then resuspend in 400 μ L of 1x Annexin V Binding Buffer for analysis, or continue to Steps 8-11 if immunophenotyping is required (optional).
- 8. **For immunophenotyping**: Resuspend cells with 50 μ L 1x Annexin V Co-staining Buffer (Item D) containing Fc Receptor blocking reagents (e.g. RayBright[®] Human or Mouse FcR Block), incubate for 5 minutes.

- 9. Add 50 μ L of Annexin V Co-staining Buffer containing your cocktail of fluorochrome-conjugated antibodies to the above cells (total volume of 100 μ L) and incubate for 25 minutes at 4°C.
- 10.Wash the cells with 200 μ L of 1x Annexin V Co-staining Buffer and spin down the cells.
- 11.Add 400 μ L of 1x Annexin V Co-staining Buffer to each well, keep samples on ice and analyze by flow Cytometry.

Representative Data:

 One million human PBMCs were stained with RayBright[®] Live R780 and RayBright[®] V450 Annexin V in 1x Annexin V Binding Buffer per the above protocol.



IV. Storage and Stability

Store kit at 4°C, protected from light for up to three months for full functionality

This product is for research use only.

