

ReadiPrep™ Dead Cell Removal Kit

Catalog number: 67300
Unit size: 1x10⁹ cells

Component	Storage	Amount (Cat No. 67300)
Component A: ReadiPrep™ Dead Cell (Annexin V) Conjugate	Refrigerated (2-8 °C)	One vial (500 µL)
Component B: Magnetic Spheres (2%)	Refrigerated (2-8 °C)	One vial (1 mL)
Component C: Isolation Buffer	Refrigerated (2-8 °C)	One bottle (10 mL)

OVERVIEW

The ReadiPrep™ Dead Cell Removal Kit provides a fast and reliable method for enriching live cell populations by removing dead cells, including both apoptotic and necrotic cells, from your samples. This kit leverages Annexin V conjugation to specifically bind dead cells, followed by magnetic separation using magnetic spheres, ensuring that live cells are efficiently isolated from dead ones.

This kit is designed for ease of use with minimal hands-on time. The provided reagents and protocol can help to quickly obtain enriched live cell populations from a variety of sample types, such as cultured cells or tissue homogenates. The resulting live cell suspension is highly pure, making it ideal for downstream applications like flow cytometry, cell culture, or molecular analysis making it ideal for studies involving cell viability, apoptosis, or cellular responses to drug treatments.

AT A GLANCE

Protocol Summary

1. Prepare cells in the Isolation Buffer (Component C).
2. Add ReadiPrep™ Dead Cell (Annexin V) Conjugate (Component A).
3. Incubate 10 min at RT.
4. Add Magnetic Spheres (Component B).
5. Incubate 10 min at RT.
6. Place the tube in a magnetic rack.
7. Wait for 2-3 min.
8. Carefully collect the supernatant containing live cells.

SAMPLE EXPERIMENTAL PROTOCOL

For processing 1x10⁷ Cells

1. Prepare 1x10⁷ cells in 500 µL Isolation Buffer (Component C).
2. Add 25 µL ReadiPrep™ Dead Cell (Annexin V) Conjugate (Component A) to the sample. Incubate for 10 min at RT.
3. Centrifuge the cells.
4. Resuspend cells in 500 µL Isolation Buffer (Component C).
5. Add 10 µL of Magnetic Spheres (Component B) to cell sample.
6. Incubate for 10 min at RT. (The dead cells will bind to the Streptavidin Magnetic Spheres).
7. Place the tube in the magnetic rack and incubate for 2 min.
8. Collect the cell suspension into a new tube (the supernatant containing live cells).

Volume of reagents are scalable for specific cell number:

Cells number	Component A	Component B	Component C
1X10 ⁷ cells	25 µL	10 µL	0.5 mL
1X10 ⁸ cells	100 µL	100 µL	2 mL
1X10 ⁹ cells	200 µL	1 mL	4 mL

EXAMPLE DATA ANALYSIS AND FIGURES

Placeholder for image details

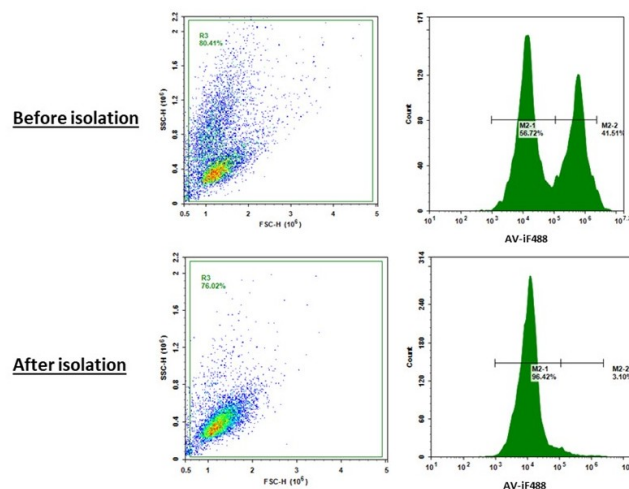


Figure 1. Starting with 72 hours cultured Jurkat cells, which has 56% live cells and 41.5% apoptosis cells (Annexin V positive), the live cell content is enriched to 94% after removal of dead cells with ReadiPrep™ Dead Cell Removal Kit (Cat#67300).

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.