

ab65782

Apoptotic Cell Enrichment Kit

Instructions for Use

For the quick and simple isolation of apoptotic cells

[View kit datasheet: www.abcam.com/ab65782](http://www.abcam.com/ab65782)

(use www.abcam.cn/ab65782 for China, or www.abcam.co.jp/ab65782 for Japan)

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1.	Overview	3
2.	Protocol Summary	4
3.	Components and Storage	5
4.	Assay Protocol	6

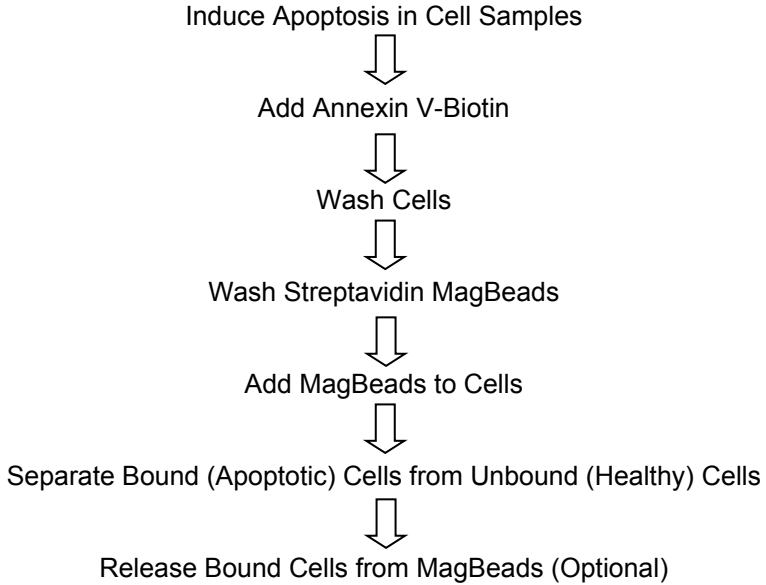
1. Overview

Abcam's Apoptotic Cell Enrichment Kit provides a simple and efficient means for isolation of apoptotic cells or removal of dead cells from cell culture or tissue preparations using annexin V magnetic beads (MagBeads).

Annexin V is a Ca^{2+} -dependent phospholipid binding protein with high affinity for phosphatidylserine (PS), which is redistributed from the inner to the outer plasma membrane leaflet in apoptotic or dead cells. Once on the cell surface, PS becomes available for binding to annexin V and any of its conjugates. Binding of annexin V-biotin to apoptotic cells followed by binding of the biotin to streptavidin-MagBeads enables separation of apoptotic cells from living cells. The apoptotic cells bound to the MagBeads adhere to the magnet, while non-apoptotic cells stay in suspension. The separated apoptotic and healthy cells can then be used in a variety of assays to study apoptotic mechanisms and pathways.

The kit has also been successfully used to remove dead cells from healthy cell culture.

2. Protocol Summary



3. Components and Storage

A. Kit Components

Item	Quantity
Annexin V-Biotin	150 µL
1X Binding Buffer	25 mL
Streptavidin MagBeads	1.5 mL
Apoptotic Cell Releasing Buffer	10 mL

* Store kit at -20°C.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Magnetic separator
- Orbital shaker

4. Assay Protocol

1. Induce apoptosis by desired method.
2. Collect $5\text{-}10 \times 10^6$ cells by centrifugation.
3. Re-suspend cells in 100 μl of 1X Binding Buffer.
4. Add 5 μl of Annexin V-Biotin, mix gently.
5. Incubate at room temperature for 5-10 min.
6. Centrifuge the cells for 2 min at 600 x g, remove the supernatant.
7. Wash the cells with 200 μl 1X Binding Buffer, repeat Step 6.
8. Re-suspend the cells in 200 μl 1X Binding Buffer.
9. Wash the Streptavidin MagBeads with 1X Binding Buffer:
 - a) Transfer 50 μl /assay of the MagBeads suspension to a new tube.
 - b) Separate the beads by a magnetic separator and remove the solvent.
 - c) Add 200 μl 1X Binding Buffer to the Beads.
 - d) Separate the Beads again and re-suspend them in 50 μl /assay 1X Binding Buffer.

- 10.** Add 50 µl of the re-suspended MagBeads to the cells.
- 11.** Rotate for 15 min at +4°C.
- 12.** Separate the MagBeads using the magnetic separator. Wait a few minutes for the separation to progress.
- 13.** Carefully transfer the unbound cells from the beads to a new tube.
- 14.** Spin the unbound cells (healthy cell population) and remove the supernatant. Keep the healthy cell population for further study.
- 15.** The magnetic bound apoptotic cells can be used directly for apoptosis assays (e.g., caspase activity assays or others) or released from the beads using the following procedure:
 - a)** Add 100 µl of the Apoptotic Cell Releasing Buffer. Gently mix and incubate for 10 min at room temp. Separate the beads using the Magnetic Separator and carefully transfer the solution (containing apoptotic cells) to a new tube.
 - b)** Repeat the releasing process again and combine the released cells together.
 - c)** Spin down the apoptotic cells and save for further study.

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “*contact us*” on www.abcam.com for the phone number for your region).

UK, EU and ROW

Email: technical@abcam.com

Tel: +44 (0)1223 696000

www.abcam.com

US, Canada and Latin America

Email: us.technical@abcam.com

Tel: 888-77-ABCAM (22226)

www.abcam.com

China and Asia Pacific

Email: hk.technical@abcam.com

Tel: 400 921 0189 / +86 21 2070 0500

www.abcam.cn

Japan

Email: technical@abcam.co.jp

Tel: +81-(0)3-6231-0940

www.abcam.co.jp