

# Apoptotic Cell Enrichment Kit ab65782

## Overview

<b>Product name</b>	Apoptotic Cell Enrichment Kit
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Direct
<b>Assay time</b>	1h 00m
<b>Product overview</b>	<p>Abcam's Apoptotic Cell Isolation 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 <math>Ca^{2+}</math>-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.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K258 Apoptotic Cell Isolation Kit. K258-30 is the same size as the 30 test size of ab65782.</p>
<b>Tested applications</b>	<b>Suitable for:</b> FM, Flow Cyt

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	30 tests
1x Binding Buffer	1 x 25ml
Annexin V-Biotin	1 x 150µl
Apoptotic Cell Releasing Buffer	1 x 10ml
Streptavidin-MagBeads	1 x 1.5ml

**Relevance** Cell isolation is a way of dissociating cells from tissues and separating the various types in order

to obtain a relatively homogeneous population of cells for analysis.

## Applications

---

### The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab65782 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
FM		Use at an assay dependent concentration.
Flow Cyt		Use at an assay dependent concentration.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors