

Apoptotic Cell Enrichment Kit ab65782

概述

产品名称	Apoptotic Cell Enrichment试剂盒
样品类型	Adherent cells, Suspension cells
检测类型	Direct
检测时间	1h 00m
产品概述	<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 Ca²⁺-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 FAQs page for tips and troubleshooting.</p>
说明	<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>
经测试应用	适用于: FM, Flow Cyt

性能

存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	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

相关性	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.

应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab65782于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
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.cn/abpromise> or contact our technical team.

Terms and conditions

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