

Annexin V- DY-634 PI Apoptosis Staining / Detection Kit ab214484

[2 References](#) [2 图像](#)

概述

产品名称	Annexin V- DY-634 PI Apoptosis Staining / Detection试剂盒
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based (quantitative)
种属反应性	与反应: Mammals, Other species
产品概述	<p>Annexin V- DY-634 PI Apoptosis Staining / Detection Kit (ab214484) contains Annexin V labeled with DY-634, which allows the identification and quantitation of apoptotic cells on a single-cell basis by flow cytometry. Simultaneous staining of cells with Annexin V – DY634 (red fluorescence) and the non-vital dye propidium iodide (PI) (orange fluorescence) allows the discrimination of intact cells (AnnexinV-DY-634 negative, PI Staining Solution negative), early apoptotic (Annexin V-DY-634 positive, PI Staining Solution negative) and late apoptotic or necrotic cells (Annexin V- DY-634 positive, PI Staining Solution positive).</p> <p>DY-634 can be used as direct replacement for APC.</p> <p>For more apoptosis assays, review the full set of Annexin V assays, or the apoptosis assay and apoptosis marker guide.</p>

说明

Apoptosis is a regulated process of cell death that occurs during embryonic development as well as maintenance of tissue homeostasis. Inappropriately regulated apoptosis is implicated in different disease states, such as neurodegeneration disease and cancer. The apoptosis program is characterized by morphologic features, including loss of plasma membrane asymmetry and attachment, condensation of the cytoplasm and nucleus, and compaction and fragmentation of the nuclear chromatin. Exposure of phosphatidylserine (PS) on the external surface of the cell membrane has been reported to occur in the early phases of apoptotic cell death, during which the cell membrane remains intact. In leukocyte apoptosis, PS on the outer surface of the cell marks the cell for recognition and phagocytosis by macrophages. The human vascular anticoagulant, annexin V, is a 35-36 kDa Ca^{2+} dependent phospholipid binding protein that has a high affinity for PS, and shows minimal binding to phosphatidylcholine and sphingomyelin. Changes in PS asymmetry, which can be analyzed by measuring annexin V binding to the cell membrane, are generally observed before morphological changes associated with apoptosis occurred and before membrane integrity is lost.

平台

Flow cytometer

性能

存放说明

Store at +4°C. Please refer to protocols.

组件	100 tests
Annexin V DY-634 Conjugate (in PBS, pH 7.4)	1 x 500µl
Propidium Iodide Staining Solution (in PBS, pH 7.4)	1 x 500µl
10X Binding Buffer Concentrate	1 x 50ml

功能

This protein is an anticoagulant protein that acts as an indirect inhibitor of the thromboplastin-specific complex, which is involved in the blood coagulation cascade.

疾病相关

Pregnancy loss, recurrent, 3

序列相似性

Belongs to the annexin family.
Contains 4 annexin repeats.

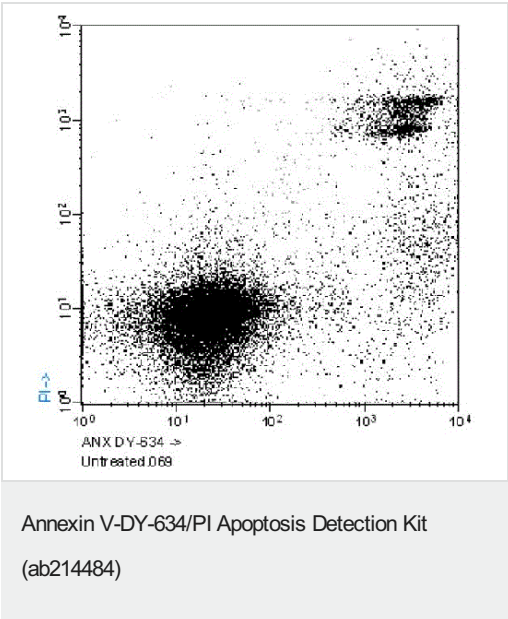
结构域

The [IL]-x-C-x-x-[DE] motif is a proposed target motif for cysteine S-nitrosylation mediated by the iNOS-S100A8/A9 transnitrosylase complex.
A pair of annexin repeats may form one binding site for calcium and phospholipid.

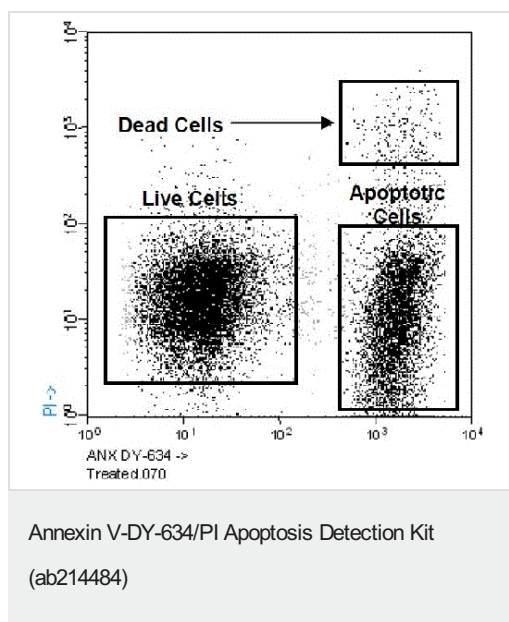
翻译后修饰

S-nitrosylation is induced by interferon-gamma and oxidatively-modified low-density lipoprotein (LDL(ox)) possibly implicating the iNOS-S100A8/9 transnitrosylase complex.

图片



Untreated Jurkat cells were stained with Annexin V-DY-634/PI Apoptosis Detection Kit (ab214484), following assay protocol.



Jurkat cells treated with 6 μ M camptothecin for four hours were stained with Annexin V-DY-634/PI Apoptosis Detection Kit (ab214484), following assay protocol. Gating showing the different populations of live, dead and apoptotic cells are shown in the figure.

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