

Annexin V-FITC Apoptosis Staining / Detection Kit ab14085

★★★★★ [3 Abreviews](#) [478 References](#) [4 图像](#)

概述

产品名称	Annexin V-FITC Apoptosis Staining / Detection试剂盒
样品类型	Adherent cells, Suspension cells
检测类型	Direct
检测时间	0h 10m
产品概述	<p>Annexin V-FITC Apoptosis Staining / Detection Kit ab14085 is used in a 10 min, one-step staining procedure to detect apoptosis by staining phosphatidylserine molecules which have translocated to the outside of the cell membrane. Analysis is by flow cytometry or fluorescence microscopy.</p> <p>The kit can differentiate apoptosis vs necrosis when performing both Annexin V-FITC and PI staining.</p> <p>The Annexin V-FITC reagent contained in the kit is also available as Annexin V-FITC reagent ab14082.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K101 Annexin V-FITC Apoptosis Kit. K101-100 is the same size as the 100 test size of ab14085.</p> <p>Soon after initiating apoptosis, cells translocate membrane phosphatidylserine molecules from the inner face of the plasma membrane to the cell surface. Phosphatidylserine on the cell surface is detected by staining with a fluorescent conjugate of Annexin V, a protein that has a high affinity for phosphatidylserine.</p> <p>For more apoptosis assays, review the full set of Annexin V assays, or the apoptosis assay and apoptosis marker guide.</p>
平台	Flow cytometer, Fluorescence microscope

性能

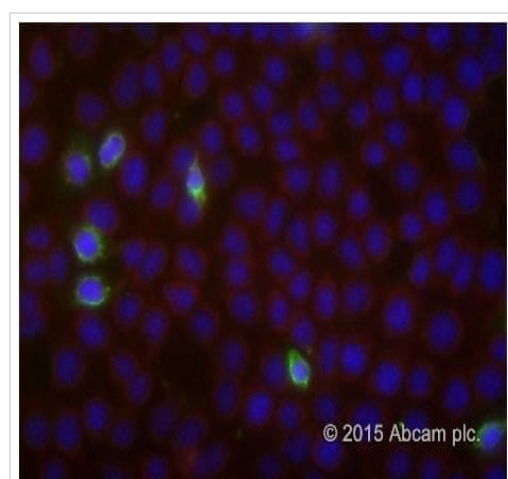
存放说明 Store at +4°C. Please refer to protocols.

组件	100 tests
Annexin V-FITC II	1 x 500µl

组件	100 tests
Binding Buffer II	1 x 50ml
Propidium Iodide II	1 x 500µl

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.

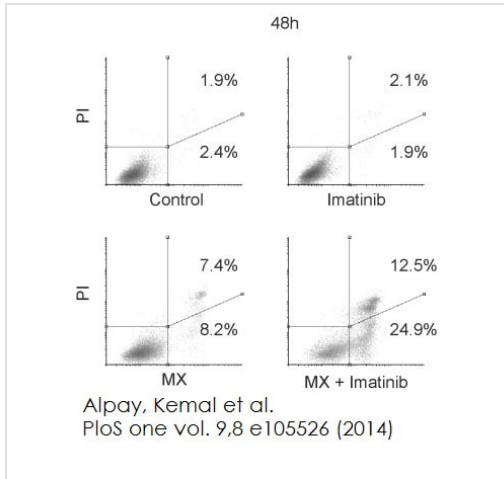
图片



Ab14085 was used to determine minor levels of apoptosis (using both the Annexin V-FITC and PI) in mouse cortical collecting duct cells (mCCDs). mCCD cells were incubated with serum free medium for 48h. The green label on the plasma membrane (Annexin V-FITC) and the absence of nuclear red (PI) staining indicates apoptosis rather than necrosis. Fluorescent microscopy was used to analyse the cells.

Apoptosis in Mouse Cortical Collecting Duct Cells

Image courtesy of an anonymous abreview

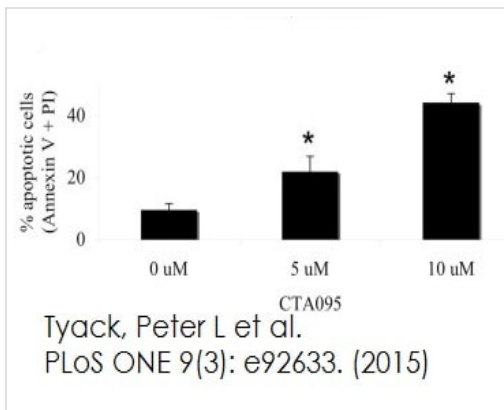


Flow cytometry analysis of treated HeLa cells for
48 hours

Alpay et al., PLoS One, 9(19), Fig 5B.
Doi:10.1371/journal.pone.0105526 Reproduced under
the Creative Commons license
<http://creativecommons.org/licenses/by/4.0/>

HeLa cells were harvested with trypsinization together with floating non-viable cells. Cells were washed with PBS and suspended in sodium citrate buffer 20 minutes prior to analysis. HeLa cells were treated with Mitoxantrone (MX) and MX +Imatinib for 48 hours. The samples were then stained with Annexin V-FITC Apoptosis Staining/Detection kit (ab14085). A FACSCalibur flow cytometer was used for cell cycle analysis.

This is a modified version of the original image

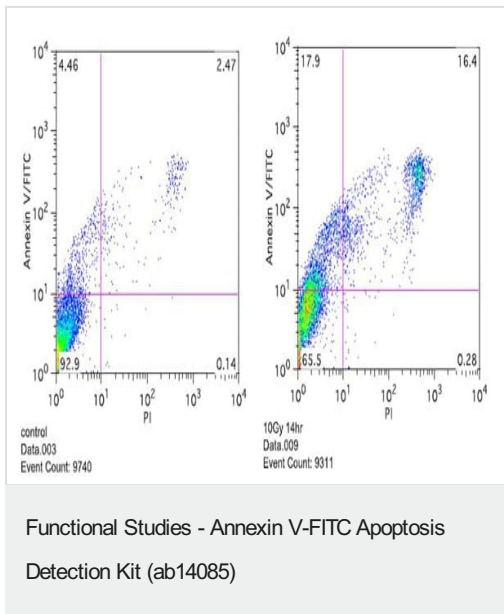


Analysis of apoptosis in prostate cancer cells
following treatment with CTA095

Guo W et al., PLoS One, 8(8). Fig7b, doi:
10.1371/journal.pone.0070910 Reproduced under the
Creative Commons license
<http://creativecommons.org/licenses/by/4.0/>

PC3 cells were seeded at 10^6 cells/ml and incubated overnight and then treated with CTA095 at various concentrations for 24hours. Apoptosis was then analyzed using Annexin-V FITC apoptosis detection kit (ab14085).

This is a modified version of the original image



Annexin V-FITC/ PI staining of AG06173 primary fibroblasts.

10^5 cells were used for analysis. Resuspended cells were incubated with Annexin V-FITC for 15 min in the dark. Propidium iodide (**ab14083**) was used as a counterstain to discriminate necrotic/ dead cells from apoptotic cells. *Left:* negative control - AG6173 untreated cells. *Right:* positive control - AG6173 cells irradiated at 10 Gy.

Image courtesy of S. Khoronenkova PhD, Gray Institute for Radiation Oncology and Biology, University of Oxford, Oxford, UK.

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