

Product datasheet

Annexin V-Cy5 Apoptosis Staining / Detection Kit ab14150

[26 References](#) [1 Image](#)

Overview

Product name	Annexin V-Cy5 Apoptosis Staining / Detection Kit
Sample type	Adherent cells, Suspension cells
Assay type	Direct
Assay time	0h 10m
Product overview	Annexin V-Cy5 Apoptosis Staining / Detection Kit ab14150 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.

Fluorescence can be detected at Ex/Em = 649/670 nm.

Analysis is by flow cytometry or fluorescence microscopy.

The Annexin V-Cy5 reagent contained in the kit is also available as Annexin V-Cy5 reagent **ab14147**.

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K103 Annexin V-Cy5 Apoptosis Kit. K103-100 is the same size as the 100 test size of ab14150.</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 <u>Annexin V assays</u>, or the <u>apoptosis assay and apoptosis marker guide</u>.</p>
--------------	--

Platform	Flow cytometer, Fluorescence microscope
-----------------	---

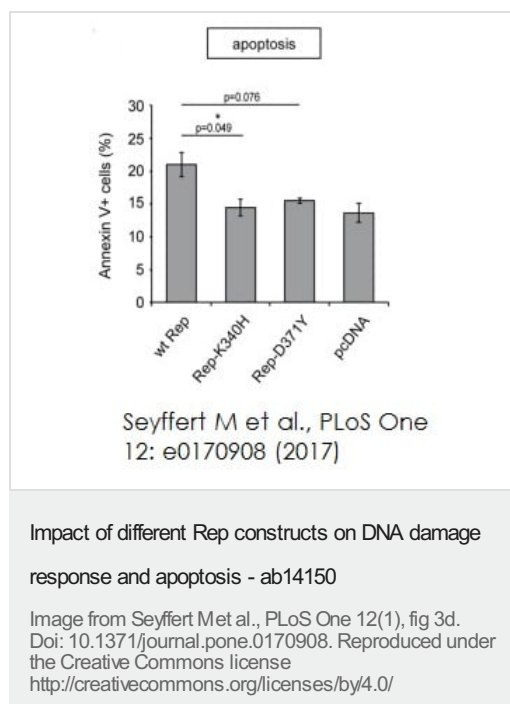
Properties

Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	100 tests
Annexin V-Cy5	1 x 500µl
Binding Buffer II	1 x 50ml

Function	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.
Involvement in disease	Pregnancy loss, recurrent, 3
Sequence similarities	Belongs to the annexin family. Contains 4 annexin repeats.
Domain	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.
Post-translational modifications	S-nitrosylation is induced by interferon-gamma and oxidatively-modified low-density lipoprotein (LDL(ox)) possibly implicating the iNOS-S100A8/9 transnitrosylase complex.

Images



Vero cells were co-transfected with different Rep-constructs together with a plasmid expressing eGFP to identify successful transfection. After three days, cells were stained using Annexin V-Cy5 Apoptosis staining/detection kit (ab14150) and analyzed by flow cytometry with filters for eGFP and Cy5.

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