

# Annexin V-EGFP Apoptosis Staining / Detection Kit ab14153

## 7 References

### 概述

产品名称	Annexin V-EGFP 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	

Annexin V-EGFP Apoptosis Staining / Detection Kit ab14153 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 with a FITC filter.

EGFP is brighter and more photo-stable than other fluorescent reagents. The kit can differentiate apoptosis vs necrosis when performing both annexin V-EGFP and PI staining.

### 说明

This product is manufactured by BioVision, an Abcam company and was previously called K104 Annexin V-EGFP Apoptosis Kit. K104-100 is the same size as the 100 test size of ab14153.

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.

For more apoptosis assays, review the full set of [Annexin V assays](#), or the [apoptosis assay and apoptosis marker guide](#).

### 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-EGFP	1 x 500µl
Binding Buffer	1 x 50ml

组件	100 tests
Propidium iodide	1 x 500µl

<b>功能</b>	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.
<b>疾病相关</b>	Pregnancy loss, recurrent, 3
<b>序列相似性</b>	Belongs to the annexin family. Contains 4 annexin repeats.
<b>结构域</b>	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.
<b>翻译后修饰</b>	S-nitrosylation is induced by interferon-gamma and oxidatively-modified low-density lipoprotein (LDL(ox)) possibly implicating the iNOS-S100A8/9 transnitrosylase complex.

**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