

# Annexin V-Biotin Apoptosis Staining / Detection Kit ab14190

## 2 References

### 概述

产品名称	Annexin V-Biotin Apoptosis Staining / Detection试剂盒
样品类型	Adherent cells, Suspension cells
检测类型	Direct
检测时间	0h 30m
产品概述	<p>Annexin V-Biotin Apoptosis Staining / Detection Kit (<b>ab1419</b>) is based on the observation that soon after initiating apoptosis, most cell types translocate the membrane phospholipid phosphatidylserine (PS) from the inner face of the plasma membrane to the cell surface. Once on the cell surface, PS can easily bind to Biotin-conjugated Annexin V, a protein that has a strong natural affinity for PS. Annexin V-Biotin can be detected in conjunction with conventional dye-staining using any streptavidine- or avidin-dye reagents, such as (strept)avidine-fluorescein, -peroxidase, -alkaline phosphatase (AP), and -beta-gal, etc.</p> <p>For more apoptosis assays, review the full set of <a href="#">Annexin V assays</a>, or the <a href="#">apoptosis assay and apoptosis marker guide</a>.</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-Biotin	1 x 500µl
Binding Buffer	1 x 50ml
Propidium iodide	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.

<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