

## Product datasheet

# Annexin V-PE-Cy5 Apoptosis Staining / Detection Kit ab14159

### Overview

<b>Product name</b>	Annexin V-PE-Cy5 Apoptosis Staining / Detection Kit
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Direct
<b>Assay time</b>	0h 10m
<b>Product overview</b>	<p>Annexin V-PE-Cy5 Apoptosis Staining / Detection Kit (ab14159) 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 be easily detected by staining with a fluorescent conjugate of Annexin V, a protein that has a strong natural affinity for PS. The one-step staining procedure takes only 10 minutes. In addition, the assay can be directly performed on live cells, without the need for fixation. Results can be analyzed by flow cytometry or fluorescence microscopy.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
<b>Platform</b>	Flow cytometer, Fluorescence microscope

### Properties

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

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

<b>Function</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>Involvement in disease</b>	Pregnancy loss, recurrent, 3
<b>Sequence similarities</b>	Belongs to the annexin family. Contains 4 annexin repeats.
<b>Domain</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.

**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.

---

**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