

## Product datasheet

# Annexin V-Cy3 Apoptosis Staining / Detection Kit ab14142

### 6 References

#### Overview

<b>Product name</b>	Annexin V-Cy3 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>	Annexin V-Cy3 Apoptosis Staining / Detection Kit ab14142 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.

The Annexin V-Cy3 reagent contained in the kit is also available as Annexin V-Cy3 reagent [ab14143](#).

<b>Notes</b>	<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 <a href="#">Annexin V assays</a>, or the <a href="#">apoptosis assay and apoptosis marker guide</a>.</p>
--------------	--

<b>Platform</b>	Flow cytometer, Fluorescence microscope
-----------------	---

#### Properties

<b>Storage instructions</b>	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	100 tests
Annexin V-Cy3	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.
<b>Post-translational modifications</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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### **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