

## Annexin V Conjugates

Annexin V fluorescent conjugates can be used to detect apoptotic cells. Annexin V is available conjugated to a wide variety of dyes.



### Product Description

Annexin V fluorescent conjugates can be used to detect apoptotic cells. The human anticoagulant Annexin V is a 35-36 kDa calcium-dependent phospholipid-binding protein with high affinity for phosphatidylserine (PS). In normal viable cells, PS is located on the inner leaflet of the cytoplasmic membrane. In apoptotic cells, however, PS is translocated from the inner to the outer leaflet of the plasma membrane, where it can be detected by fluorescently labeled Annexin V.

- Detect apoptotic cells with fast, reliable detection of phosphatidylserine externalization
- Conjugates with wide choice of CF® Dye colors, plus biotin, R-PE, APC, and other fluorescent dyes
- Biotium's CF® Dyes have superior brightness, photostability, and are water-soluble
- Can be used for fluorescence microscopy, live cell imaging, and flow cytometry

CF® Dye, biotin, and FITC conjugates of Annexin V are provided as 50 ug/mL solutions in 10 mM Tris, 1 mM EDTA, 30 mM NaCl, 1 mg/mL bovine serum albumin, 0.1% sodium azide, pH 7.5. R-PE and APC Annexin V conjugates are supplied in PBS, 5 mg/mL BSA, 0.05% sodium azide at 5 uL per test.

### Detect Apoptosis

We also offer [Annexin V Conjugates, Azide-Free and Lyophilized](#) in colors ranging from UV to far-red for live cell staining, and [Near-IR CF® Dye Annexin V](#) (also lyophilized and preservative-free), compatible with small animal NIR *in vivo* imaging. Annexin V is also available with other probes in our [Apoptotic and Necrotic Staining Kits](#). See our full selection of [Cell Viability and Apoptosis Assays](#).

### Superior CF® Dyes

Biotium's next-generation CF® Dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about CF® Dyes.

**Note:** Conjugates of blue-fluorescent dyes like CF®350, CF®405S and CF®405M are not recommended for detecting low abundance targets and may be challenging to use in tissue specimens. Blue dyes have lower fluorescence and photostability, and cells and tissue have high autofluorescence in blue wavelengths, resulting in lower signal to noise compared to other colors.

### Product attributes

<b>Storage Conditions</b>	Store at 2 to 8 °C, Do not freeze, Protect fluorescent conjugates from light
<b>Apoptosis/viability marker</b>	Phosphatidylserine/Annexin V
<b>For live or fixed cells</b>	For live/intact cells
<b>Assay type/options</b>	Endpoint assay, Homogeneous assay, No-wash staining, Real-time imaging, Short term staining (<24h)
<b>Detection method/readout</b>	Fluorescence microscopy, Live cell imaging, Flow cytometry
<b>Format</b>	Fluorescent dye & biotin conjugates: Supplied in 10 mM Tris pH 7.5, 1 mM EDTA, 30 mM NaCl, 1 mg/mL BSA, 0.1% sodium azide R-PE & APC conjugates: Supplied in PBS, 5 mg/mL BSA, 0.05% sodium azide
<b>Colors</b>	Blue, Green, Orange, Red, Far-red, Near-infrared
<b>Fixation options</b>	Fix after staining (formaldehyde)
<b>Product origin</b>	Annexin V (human); recombinant, produced in E. coli, Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

## Annexin V Conjugates

Product	Ex/Em	Concentration	Size	Catalog No.
CF@350 Annexin V	347/448 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29012</a>
CF@405M Annexin V	408/452 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29009</a>
CF@450 Annexin V	450/538 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29083</a>
CF@488A Annexin V	490/515 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29005</a>
CF@555 Annexin V	555/565 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29004</a>
CF@568 Annexin V	562/583 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29010</a>
CF@583R Annexin V	586/609 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29085</a>
CF@594 Annexin V	593/614 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29011</a>
CF@633 Annexin V	630/650 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29008</a>
CF@640R Annexin V	642/662 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29014</a>
CF@647 Annexin V	650/665 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29003</a>
CF@660R Annexin V	663/682 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29069</a>
CF@740 Annexin V	742/767 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29126</a>
R-PE Annexin V	496, 546, 565/578 nm	5 uL/test	100 uL	<a href="#">Purchase 29045-100uL</a>
500 uL	<a href="#">Purchase 29045-500uL</a>			
APC Annexin V	650/660 nm	5 uL/test	100 uL	<a href="#">Purchase 29057-100uL</a>
500 uL	<a href="#">Purchase 29057-500uL</a>			
FITC Annexin V	496/516 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29001</a>
Biotin Annexin V	N/A	50 ug/mL	0.5 mL	<a href="#">Purchase 29013</a>
5X Annexin V Binding Buffer	N/A	5X	15 mL	<a href="#">Purchase 99902</a>

CF is a registered trademark of Biotium, Inc. Alexa Fluor, and DyLight are registered trademarks of Thermo Fisher Scientific.

## References

Download a list of curated [CF@ Dye Annexin V references](#).

This datasheet was generated on June 5, 2026 at 02:54:04 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/annexin-v-conjugates/>