



Glowing products for science

## Annexin V Conjugates, Azide-Free, Lyophilized

Preservative-free Annexin V conjugates are compatible with real-time staining of apoptotic cells for live cell imaging, fluorescence microscopy, or flow cytometry. Near-IR Annexin V fluorescent conjugates are suitable for *in vivo* imaging.



### Product Description

CF® Dye Annexin V Conjugates can be used to stain the surface of 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.

- Azide-free for no-wash, real-time live cell imaging in culture medium
- Near-IR CF® Dyes are compatible with small animal *in vivo* imaging
- Fast & simple detection of phosphatidylserine on apoptotic cells
- Choice of 19 CF® Dye colors with superior brightness & photostability
- For real-time live cell imaging, fluorescence microscopy, or flow cytometry

### Preservative Free for Real-Time, Live Cell Imaging

Annexin V conjugates typically are supplied as stock solutions with azide as a preservative for end-point staining assays in Annexin V binding buffer. Our azide-free CF® Dye Annexin V Conjugates are supplied as lyophilized solids with no azide or other preservatives that might be incompatible with live cell or *in vivo* imaging. After reconstitution in buffer, the conjugates can be added to cell culture medium for no-wash, real-time live cell imaging. Our [Mini Syringe Filters](#) are convenient for small volume sterile filtration of azide-free Annexin V stock solutions or other aqueous solutions for use in cell culture.

See Annexin V staining in real time:

### 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 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.

### More Apoptosis Assays

We also offer [Annexin V Conjugate](#) solutions (with azide) with a large selection of CF® Dyes, biotin, R-PE, APC, and other labels. In addition, we offer [Apoptotic and Necrotic Staining Kits](#) containing Annexin V and other probes. See our full selection of [Cell Viability and Apoptosis Assays](#).

Call us : [800-304-5357](tel:800-304-5357)

### Product attributes

|                                   |  |
|-----------------------------------|--|
| <b>Apoptosis/viability marker</b> | Phosphatidylserine/Annexin V   |
| <b>For live or fixed cells</b>    | For live/intact cells  |
| <b>Detection method/readout</b>   | Fluorescence microscopy, Live cell imaging, Flow cytometry, Near-IR imager, <i>In vivo</i> near-IR imaging |
| <b>Assay type/options</b>         | Endpoint assay, Homogeneous assay, Long term staining (24-72h), No-wash staining, Real-time imaging        |
| <b>Fixation options</b>           | Fix after staining (formaldehyde)  |
| <b>Colors</b>                     | Blue, Green, Red, Far-red, Near-infrared   |
| <b>Product origin</b>             | Annexin V (human); recombinant, produced in <i>E. coli</i>   |

CF® Dye Annexin V Conjugates, Azide-Free, Lyophilized

| Conjugation             | Ex/Em      | Size  | Catalog No.                | Dye Features                     |
|-------------------------|------------|-------|----------------------------|----------------------------------|
| <a href="#">CF@350</a>  | 347/448 nm | 5 ug  | <a href="#">29012R-5ug</a> | <a href="#">CF@350 Features</a>  |
| <a href="#">CF@405M</a> | 408/452 nm | 5 ug  | <a href="#">29009R-5ug</a> | <a href="#">CF@405M Features</a> |
| <a href="#">CF@450</a>  | 450/538 nm | 5 ug  | <a href="#">29083R-5ug</a> | <a href="#">CF@450 Features</a>  |
| <a href="#">CF@488A</a> | 490/515 nm | 5 ug  | <a href="#">29005R-5ug</a> | <a href="#">CF@488A Features</a> |
| <a href="#">CF@555</a>  | 555/565 nm | 5 ug  | <a href="#">29004R-5ug</a> | <a href="#">CF@555 Features</a>  |
| <a href="#">CF@568</a>  | 562/583 nm | 5 ug  | <a href="#">29010R-5ug</a> | <a href="#">CF@568 Features</a>  |
| <a href="#">CF@583R</a> | 586/609 nm | 5 ug  | <a href="#">29085R-5ug</a> | <a href="#">CF@583R Features</a> |
| <a href="#">CF@594</a>  | 593/614 nm | 5 ug  | <a href="#">29011R-5ug</a> | <a href="#">CF@594 Features</a>  |
| <a href="#">CF@633</a>  | 630/650 nm | 5 ug  | <a href="#">29008R-5ug</a> | <a href="#">CF@633 Features</a>  |
| <a href="#">CF@640R</a> | 642/662 nm | 5 ug  | <a href="#">29014R-5ug</a> | <a href="#">CF@640R Features</a> |
| <a href="#">CF@647</a>  | 650/665 nm | 5 ug  | <a href="#">29003R-5ug</a> | <a href="#">CF@647 Features</a>  |
| <a href="#">CF@660R</a> | 663/682 nm | 5 ug  | <a href="#">29069R-5ug</a> | <a href="#">CF@660R Features</a> |
| <a href="#">CF@680</a>  | 681/698 nm | 25 ug | <a href="#">29007</a>      | <a href="#">CF@680 Features</a>  |
| <a href="#">CF@680R</a> | 680/701 nm | 25 ug | <a href="#">29070</a>      | <a href="#">CF@680R Features</a> |
| <a href="#">CF@700</a>  | 696/721 nm | 25 ug | <a href="#">29082</a>      | <a href="#">CF@700 Features</a>  |
| <a href="#">CF@750</a>  | 755/777 nm | 25 ug | <a href="#">29006</a>      | <a href="#">CF@750 Features</a>  |
| <a href="#">CF@770</a>  | 770/797 nm | 25 ug | <a href="#">29046</a>      | <a href="#">CF@770 Features</a>  |
| <a href="#">CF@790</a>  | 784/806 nm | 25 ug | <a href="#">29047</a>      | <a href="#">CF@790 Features</a>  |
| <a href="#">CF@800</a>  | 797/816 nm | 25 ug | <a href="#">29078</a>      | <a href="#">CF@800 Features</a>  |

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

This datasheet was generated on December 6, 2025 at 09:40:35 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/annexin-v-conjugates-azide-free-lyophilized/>