

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](#).

Product attributes

Apoptosis/viability marker	Phosphatidylserine/Annexin V
For live or fixed cells	For live/intact cells
Detection method/readout	Fluorescence microscopy, Live cell imaging, Flow cytometry, Near-IR imager, <i>In vivo</i> near-IR imaging
Assay type/options	Endpoint assay, Homogeneous assay, Long term staining (24-72h), No-wash staining, Real-time imaging
Fixation options	Fix after staining (formaldehyde)
Colors	Blue, Green, Red, Far-red, Near-infrared
Product origin	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
CF@350	347/448 nm	5 ug	29012R-5ug	CF@350 Features
CF@405M	408/452 nm	5 ug	29009R-5ug	CF@405M Features
CF@450	450/538 nm	5 ug	29083R-5ug	CF@450 Features
CF@488A	490/515 nm	5 ug	29005R-5ug	CF@488A Features
CF@555	555/565 nm	5 ug	29004R-5ug	CF@555 Features
CF@568	562/583 nm	5 ug	29010R-5ug	CF@568 Features
CF@583R	586/609 nm	5 ug	29085R-5ug	CF@583R Features
CF@594	593/614 nm	5 ug	29011R-5ug	CF@594 Features
CF@633	630/650 nm	5 ug	29008R-5ug	CF@633 Features
CF@640R	642/662 nm	5 ug	29014R-5ug	CF@640R Features
CF@647	650/665 nm	5 ug	29003R-5ug	CF@647 Features
CF@660R	663/682 nm	5 ug	29069R-5ug	CF@660R Features
CF@680	681/698 nm	25 ug	29007	CF@680 Features
CF@680R	680/701 nm	25 ug	29070	CF@680R Features
CF@700	696/721 nm	25 ug	29082	CF@700 Features
CF@750	755/777 nm	25 ug	29006	CF@750 Features
CF@770	770/797 nm	25 ug	29046	CF@770 Features
CF@790	784/806 nm	25 ug	29047	CF@790 Features
CF@800	797/816 nm	25 ug	29078	CF@800 Features

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

This datasheet was generated on June 4, 2026 at 05:34:29 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/annexin-v-conjugates-azide-free-lyophilized/>