



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

Annexin V CF® Dye 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.



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
- Fast & simple detection of phosphatidylserine on apoptotic cells
- Choice of 11 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 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, 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.

More Apoptosis Assays

We also offer [Near-IR CF® Dye Annexin V Conjugates](#), preservative-free and lyophilized, compatible with small animal in vivo imaging. [Annexin V Conjugate](#) solutions (with azide) also are available with a large selection of CF® Dyes, biotin, R-PE, APC, and other labels. 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](#).

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
Assay type/options	Endpoint assay, Homogeneous assay, Long term staining (24-72h), No-wash staining, Real-time imaging
Colors	Blue, Green, Red, Far-red
Fixation options	Fix after staining (formaldehyde)
Product origin	Annexin V (human); recombinant, produced in E. coli

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	5 ug	29012-5ug	CF@350 Features
CF@405M	408/452 nm	5 ug	29009-5ug	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 is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

References

Download a list of [CF® dye Annexin V references](#).

This datasheet was generated on November 22, 2024 at 11:59:42 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-annexin-v-conjugates-azide-free/>