Glowing products for science

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.

ulotum

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.

Call us: 800-304-5357 Email: btinfo@biotium.com

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
<u>CF®450</u>	450/538 nm	5 ug	29083R-5ug	CF®450 Features
<u>CF®488A</u>	490/515 nm	5 ug	29005R-5ug	CF®488A Features
<u>CF®555</u>	555/565 nm	5 ug	29004R-5ug	CF®555 Features
<u>CF®568</u>	562/583 nm	5 ug	29010R-5ug	CF®568 Features
<u>CF®583R</u>	586/609 nm	5 ug	29085R-5ug	CF®583R Features
<u>CF®594</u>	593/614 nm	5 ug	29011R-5ug	CF®594 Features
<u>CF®633</u>	630/650 nm	5 ug	29008R-5ug	CF®633 Features
<u>CF®640R</u>	642/662 nm	5 ug	29014R-5ug	CF®640R Features
CF®647	650/665 nm	5 ug	29003R-5ug	CF®647 Features
<u>CF®660R</u>	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 4, 2025 at 06:01:38 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cf-dye-annexin-v-conjugates-azide-free/