

Annexin V Near IR CF® Dye Conjugates

Near-IR Annexin V fluorescent conjugates can be used to detect apoptotic cells and are suitable for *in vivo* imaging.



Product Description

Annexin V is a 35-36 kDa protein that has a high affinity for phosphatidylserine (PS). In normal viable cells, PS is located on the cytoplasmic surface of the cell membrane. In apoptotic cells, however, PS is translocated from the inner to the outer leaflet of the plasmic membrane, making itself available for Annexin V binding. Near-IR CF® Dye Annexin V does not contain azide or other potentially toxic materials, and thus is suitable for *in vivo* imaging.

- Detect apoptotic cells with fast, reliable detection of phosphatidylserine externalization
- No azide or other potentially toxic materials or preservatives, suitable for *in vivo* imaging
- Choice of 7 near-IR fluorescent CF® Dye colors
- Biotium's CF® Dyes are far more photostable and brighter than other similar near-IR dyes
- Can be used for fluorescence microscopy, live cell imaging, and flow cytometry

Superior CF® Dyes

CF® Dyes are far more photostable and brighter than other similar near-IR dyes, such as Alexa Fluor® 700, Alexa Fluor® 750, Cy®7 and Iridye® 800. Learn more about Near-IR CF® Dyes in the [Near-IR Dye Flyer](#). Learn more about [CF® Dyes](#).

Detect Apoptosis with Annexin V

We also offer [Annexin V Conjugates, Azide-Free, and Lyophilized](#) in colors ranging from UV to far-red for live cell staining. [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](#).

Annexin V Near IR CF® Dye Conjugates

Conjugation	Ex/Em	Size	Catalog No.	Dye 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	695/720 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, 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 5, 2025 at 06:16:22 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/annexin-v-near-ir-cf-dye-conjugates/>

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, In vivo 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	Near-infrared
Product origin	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.