

## 5X Annexin V Binding Buffer

Concentrated buffer used for Annexin V staining of cells to monitor apoptosis.



### Product attributes

Storage Conditions

Store at 2 to 8 °C

## Product Description

5X Annexin V Binding Buffer is used to promote optimal Annexin V staining of apoptotic cells.

- Buffer optimized for Annexin V staining of apoptotic cells
- Provided as a 5X concentrate
- Dilute to 1X in water for use during and after Annexin V staining

## Detect Apoptosis with Annexin V

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.

Biotium offers Annexin V conjugated to a [wide range of CF® Dyes](#) and other fluorescent dyes, including our unique range of [near-infrared CF® Dyes](#). We also provide [azide-free Annexin V conjugates](#) for live cell staining.

## Annexin V conjugates

Product	Ex/Em	Concentration	Size	Catalog No.
CF@350 Annexin V	347/448 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29012</a>
CF@405M Annexin V	408/452 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29009</a>
CF@450 Annexin V	450/538 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29083</a>
CF@488A Annexin V	490/515 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29005</a>
CF@555 Annexin V	555/565 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29004</a>
CF@568 Annexin V	562/583 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29010</a>
CF@583R Annexin V	586/609 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29085</a>
CF@594 Annexin V	593/614 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29011</a>
CF@633 Annexin V	630/650 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29008</a>
CF@640R Annexin V	642/662 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29014</a>
CF@647 Annexin V	650/665 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29003</a>
CF@660R Annexin V	663/682 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29069</a>
CF@740 Annexin V	742/767 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29126</a>
R-PE Annexin V	496, 546, 565/578 nm	5 uL/test	100 uL	<a href="#">Purchase 29045-100uL</a>
500 uL	<a href="#">Purchase 29045-500uL</a>			
APC Annexin V	650/660 nm	5 uL/test	100 uL	<a href="#">Purchase 29057-100uL</a>
500 uL	<a href="#">Purchase 29057-500uL</a>			
FITC Annexin V	496/516 nm	50 ug/mL	0.5 mL	<a href="#">Purchase 29001</a>
Biotin Annexin V	N/A	50 ug/mL	0.5 mL	<a href="#">Purchase 29013</a>
5X Annexin V Binding Buffer	N/A	5X	15 mL	<a href="#">Purchase 99902</a>

## References

See a list of CF® Dye Annexin V references in the Supporting Documents tab.

This datasheet was generated on June 5, 2026 at 08:14:00 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5x-annexin-v-binding-buffer/>