



Glowing products for science

## CF®488A Annexin V and 7-AAD Apoptosis Kit

This kit contains CF®488A Annexin V to stain apoptotic cells, and 7-AAD to stain necrotic cells for flow cytometry.



### Product Description

This kit contains CF®488A Annexin V for staining apoptotic cells green, and 7-AAD for staining necrotic cells with far-red fluorescence, for detection by flow cytometry.

In apoptotic cells, the phospholipid phosphatidylserine (PS) is translocated from the inner to the outer surface of the plasma membrane, which targets the dying cells for phagocytosis. The human anticoagulant, Annexin V, is a 35 kDa  $\text{Ca}^{2+}$ -dependent phospholipid binding protein with a high affinity for PS. Annexin V labeled with CF®488A labels apoptotic cells with green fluorescence by binding to PS exposed on the outer leaflet. Our CF®488A dye is superior to fluorescein/FITC because it is not affected by pH and has far better photostability.

7-AAD is a membrane-impermeant nucleic acid dye that is excluded by live cells, but stains necrotic cells and late apoptotic cells. 7-AAD (Abs/Em with DNA = 543/655 nm) emits in the far-red spectrum and can be well separated from CF®488A and other green emitting dyes using the 488 nm line of a flow cytometer.

See our full selection of [Cell Viability & Apoptosis Assays](#).

CF dye technology is covered by US and international patents. Cy Dye is a registered trademark of Cytiva.

### References

Download a list of curated [CF® Dye references](#).

This datasheet was generated on January 8, 2026 at 01:10:11 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cf488a-annexin-v-and-7-aad-apoptosis-kit/>

Call us : [800-304-5357](tel:800-304-5357)

### Product attributes

<b>Apoptosis/viability marker</b>	Phosphatidylserine/Annexin V, Dead cell stain, Apoptosis/necrosis assay
<b>For live or fixed cells</b>	For live/intact cells
<b>Detection method/readout</b>	Flow cytometry
<b>Assay type/options</b>	Endpoint assay
<b>Colors</b>	Green/Far-red
<b>Product origin</b>	Annexin V (human); recombinant, produced in E. coli
<b>Storage Conditions</b>	Store at 2 to 8 °C, Do not freeze, Protect from light