



Glowing products for science

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

Recombinant Annexin V (Lyophilized)

Recombinant human Annexin V expressed from E. coli as a single, non-glycosylated polypeptide chain. The protein is available with no tag or a C-terminal 6X His tag.



Product attributes

Apoptosis/viability marker	Phosphatidylserine/Annexin V
Molecular weight	36.7 kDa (No Tag), 37.6 kDa (His Tag)
Product origin	Annexin V (human); recombinant, produced in E. coli
Storage Conditions	Store at -10 to -35 °C

Product Description

Recombinant human Annexin V expressed from E. coli as a single, non-glycosylated polypeptide chain. The protein is available with no tag or a C-terminal 6X His tag.

Annexin V is a member of the annexin family of calcium-dependent phospholipid binding proteins with high affinity for phosphatidylserine (PS). In viable cells, PS is predominantly located on the inner leaflet of the plasma membrane. During apoptosis, PS translocates from the inner leaflet to the outer leaflet. Fluorescently labeled Annexin V can bind to PS located on the outer leaflet and therefore is commonly used as a marker of apoptosis. Annexin V has also been identified as an anticoagulant protein that is involved in the blood coagulation cascade through indirect inhibition of the thromboplastin-specific complex.

We also offer [Annexin V Conjugates](#) and [Annexin V Conjugates, Azide-Free and Lyophilized](#) in colors ranging from UV to far-red for live cell staining. Also view our [Annexin V Near-IR CF® Dye Conjugates](#) (also lyophilized and preservative-free), compatible with small animal NIR *in vivo* imaging. 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](#).

This datasheet was generated on January 2, 2026 at 09:17:13 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/annexin-v-his-tag/>