

# PRODUCT INFORMATION



## Annexin V (human, recombinant)

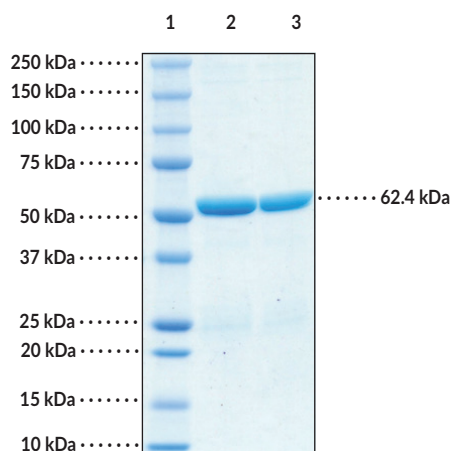
Item No. 20594

### Overview and Properties

**Synonyms:** Anchorin CII, Annexin A5, Calphobindin I, Endonexin II, Lipocortin V, PP4  
**Source:** Active recombinant N-terminal GST-tagged annexin V expressed in *E. coli*  
**Amino Acids:** 2-320 (full length)  
**Uniprot No.:** P08758  
**Molecular Weight:** 62.4 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** *batch specific*  
**Supplied in:** PBS, pH 7.4, with 10% glycerol  
**Protein**  
**Concentration:** *batch specific* mg/ml  
**Activity:** Suitable for cell-based assays

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: Annexin V (4 µg)  
Lane 3: Annexin V (2 µg)

Representative gel image shown; actual purity may vary between each batch.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/16/2020

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Annexin V is a soluble membrane-binding protein encoded by the ANXA5 gene in humans that binds to negatively charged phospholipids, such as phosphatidylserine (PS), in a calcium-dependent manner.<sup>1-3</sup> It is expressed in a wide variety of tissues but is detected primarily in placenta and vascular endothelium.<sup>4</sup> Annexin V is comprised of a variable N-terminal domain and four highly conserved annexin repeats that form the membrane-binding core.<sup>1,5</sup> Upon membrane binding, annexin V forms a two-dimensional lattice that is stabilized by protein-protein interactions.<sup>2,5</sup> In activated platelets and apoptotic cells, plasma membrane asymmetry, but not integrity, is disrupted and PS becomes exposed at the cell surface, following which annexin V binds to the surface-exposed PS and inhibits its procoagulant and pro-inflammatory activities.<sup>2,4,6,7</sup> Because of its affinity for PS in apoptotic cell membranes, fluorescent and affinity-labeled versions of annexin V have been used in affinity assays for the detection and quantification of apoptotic cells.<sup>7</sup> Annexin V levels are elevated in tumor tissue from patients with a variety of cancers, including follicular thyroid adenoma, papillary thyroid carcinoma, cervical squamous cell carcinoma, and colorectal cancer.<sup>5</sup> Cayman's Annexin V (human, recombinant) protein can be used for Western blot, ELISA, and cell-based assays.

## References

---

1. Bouter, A., Carmelle, R., Gounou, C., *et al.* Review: Annexin-A5 and cell membrane repair. *Placenta* **36(Suppl. 1)**, S43-S49 (2015).
2. Reutelingsperger, C.P. and van Heerde, W.L. Annexin V, the regulator of phosphatidylserine-catalyzed inflammation and coagulation during apoptosis. *Cell. Mol. Life Sci.* **53(6)**, 527-532 (1997).
3. Bodganova, N., Horst, J., Chlystun, M., *et al.* A common haplotype of the annexin A5 (ANXA5) gene promoter is associated with recurrent pregnancy loss. *Hum. Mol. Genet.* **16(5)**, 573-578 (2007).
4. Esposito, G., Tamby, M.C., Chanseaud, Y., *et al.* Anti-annexin V antibodies: Are they prothrombotic? *Autoimmun. Rev.* **4(1)**, 55-60 (2005).
5. Peng, B., Guo, C., Guan, H., *et al.* Annexin A5 as a potential marker in tumors. *Clin. Chim. Acta* **427**, 42-48 (2014).
6. Andree, H.A., Stuart, M.C., Hermens, W.T., *et al.* Clustering of lipid-bound annexin V may explain its anticoagulant effect. *J. Biol. Chem.* **267(25)**, 17907-17912 (1992).
7. van Engeland, M., Nieland, L.J.W., Ramaekers, F.C.S., *et al.* Annexin V-affinity assay: A review on an apoptosis detection system based on phosphatidylserine exposure. *Cytometry* **31(1)**, 1-9 (1998).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM