

## Annexin V VHH Antibody

Z-N002

Support: echelon@echelon-inc.com



### **Description:**

Recombinant anti-human Annexin V VHH Antibody was raised in llama and purified from E. coli as a VHH single domain antibody with a COOH-terminal HA epitope tag.

### **Applications:**

Pull-down (PD), Immunoprecipitation (IP), and Enzyme-Linked Immunoassay (ELISA).

Other in vitro, in vivo, cellular, & tissue applications are possible, but have not been verified by Echelon Biosciences.

### **Properties:**

Form – liquid

Storage instructions – Store at -20 °C upon arrival. Aliquot and store at -80 °C if longer storage is necessary. Avoid repeated freeze/thaw cycles.

Storage buffer – 20 mM Tris-HCl pH 8.0, 150 mM NaCl, 1mM DTT, 60 % glycerol

Concentration – 1.0 mg/mL

Purity – >90%

Immunogen – purified full-length human recombinant protein

Clonality – monoclonal

Isotype – VHH single domain antibody

Tag – C-Terminal HA

### **Specificity:**

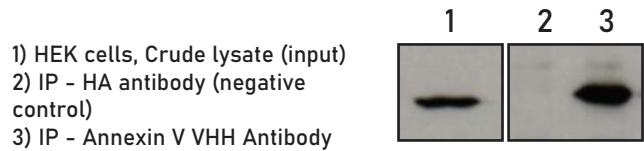
Z-N002 reacts with human Annexin V. Other species have not been tested.

### **Background:**

Annexin V, aka Annexin A5, is a 36 kDa protein that binds to the negatively charged phospholipid phosphatidylserine, a marker of apoptosis when it is exposed on the plasma membrane outer leaflet. Annexin V is often used as an indicator of apoptosis in flow cytometry experiments.

### **Data: Immunoprecipitation**

Blot:  
Anti-Annexin V antibody



HEK293T cell lysate was incubated with control antibody or Annexin V VHH Antibody followed by SDS-PAGE and WB with a mouse monoclonal Annexin V antibody.

### **References:**

1. Steels, A. et al. (2018). Intracellular displacement of p53 using transactivation domain (p53 TAD) specific nanobodies. *mAbs*; 10(7): 1045–1059.

### **Related Products:**

Products	Catalog Number
Assays and Reagents	
PI(4,5)P2 Mass Assay	K-4500
Phosphatidic Acid Beads	P-B0PA
Nanobodies	
Pan-Akt VHH Antibody	Z-N001
p53 DBD VHH Antibody	Z-N006
Tensin 4 VHH Antibody	Z-N010

Technical Data Sheet Rev. 1a, 08-31-21 - For research use only. Not intended or approved for diagnostic or therapeutic use.