

---

## Annexin V Tag Antibody

(Catalog No. A26006)

---

### Description

Mouse Anti Human Annexin V specifically recognizes human Annexin V, a phospholipid binding protein which binds with high affinity and selectivity to phosphatidylserine (PS), in the presence of calcium. The translocation of PS from the cytosolic leaflet to the outer leaflet of the plasma membrane, due to a loss in plasma membrane asymmetry during apoptosis, acts as an in vivo signal for the removal of the dying cell by phagocytosis. Detection of PS using Annexin V therefore provides a reliable diagnostic tool for the indication of early programmed cell death. Annexin V has also been identified as an anticoagulant protein in the blood coagulation cascade, by acting as an inhibitor of prothrombin activation. The presence of antibodies to Annexin V are associated with systemic lupus erythematosus (SLE), recurrent spontaneous abortions, and systemic sclerosis (SSc). Clone VAA-33 has been reported to work in western blotting applications.

### Antibody Type

Monoclonal Antibody

### Purification

Purified IgG - liquid

### Immunogen

Recombinant Annexin V

### Isotype

IgG2a

### Size

100 µg

### Formulation

Phosphate buffered saline. 0.09% Sodium Azide (NaN<sub>3</sub>)

### Specificity

Human

### Storage

Store at +4°C or at -20°C if preferred. Storage in frost-free freezers is not recommended. This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody.

### Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

### Applications & Suggested Dilutions

ELISA; Flow Cytometry: (Use 10ul of the suggested working dilution to label 1x10<sup>6</sup> cells in 100ul. Method sheets are available on request).

### Ordering Information:

Products	Size	Cat. No.
Annexin V Tag Antibody	100 µg	A26006-1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*