

Annexin V

Catalog No.	Amount	Storage	Stability
A029A	500 µg	<ul style="list-style-type: none"> • 2-6 °C • Reconstituted Annexin V should be stored in working aliquots at -20°C 	The product is stable for 1 year when stored as directed.
A029B	1 mg		

Product Description

Annexin V is a family of calcium-dependent phospholipid-binding proteins, which bind to phosphatidylserine (PS) to identify apoptotic cells. In healthy cells, PS is predominantly located along the cytosolic side of the plasma membrane. Upon initiation of apoptosis, PS loses its asymmetric distribution in the phospholipid bilayer and translocates to the extracellular membrane, which is detectable with fluorescently labeled Annexin V. In early stages of apoptosis, the plasma membrane excludes viability dyes such as propidium iodide (PI) and 7-AAD, therefore cells which display only Annexin V staining (PI/7-AAD negative) are in early stages of apoptosis. During late-stage apoptosis, loss of cell membrane integrity allows Annexin V binding to cytosolic PS, as well as cell uptake of PI and 7-AAD. Annexin V staining, paired with 7-AAD or PI is widely used to identify apoptotic stages by flow cytometry.

Source: E. coli
Molecular Weight: 35.8 kDa
Form: Lyophilized
Purity: Annexin V is >90% pure as demonstrated by SDS-gel electrophoresis.
Reconstitution: We recommend a quick spin followed by reconstitution in 1 ml distilled water (incubate at 30°C for 5 minutes). Reconstituted solution contains 20 mM phosphate buffer (pH 7.0).