

Instruction Manual

Catalog Number	PK-AB577-3357
Description	The annexin family of calcium-binding proteins is composed of at least ten mammalian genes. It is characterized by a conserved core domain which binds to phospholipids in a Calcium dependent manner and a unique amino terminal region which may confer binding specificity. Annexin family members have been implicated as regulators of such diverse processes as ion flux, endocytosis and exocytosis, and cellular adhesion. Annexin V is ubiquitously expressed at high levels in tissues and cells grown in culture.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Recombinant human Annexin V (ID BV-201)
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody detects Annexin V from samples of human, mouse, and rat origin.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-Annexin V polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The antibody can be used for Western blotting (0.5-4 µg/ml). Based on researcher's feedback, it can also be used in immunoprecipitation, and immunohistochemistry. However, the optimal conditions should be determined individually. Jurkat cell lysate (Cat.# PK-AB718-1205), mouse small intestine lysate (Cat.# PK-AB718-1408) and rat kidney tissue lysate (Cat.# PK-AB718-1465) can be used as positive controls. Blocking peptide is available separately (Cat.# PK-AB577-3357P).

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.