

Instruction Manual

Catalog Number	PK-CA577-1005-1000
Description	During apoptosis, phosphatidylserine (PS) is translocated from the cytoplasmic face of the plasma membrane to the cell surface. Annexin V has a strong, Ca ²⁺ -dependent affinity for PS and therefore is used as a probe for detecting apoptosis. The unlabeled Annexin V reagent can be conjugated to any tag of choice for detection of apoptosis. The reagent can also be used in competition binding experiments with tagged annexin V conjugates. Annexin V-FITC, -Cy3, -Cy5, -EGFP, -PE, -Biotin Labels are also available.
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	NA
Source	<i>E. coli</i>
Purity	>98% as determined by SDS-PAGE.
Appearance / Formulation / Solubility	Lyophilized. Reconstitute in water or PBS to a final concentration of 1-5 mg/ml.
Storage & Stability	Store at 4°C (short-term) or -20°C (long-term). Stable for 1 year under proper conditions.
Applications	see Description
References	NA
Caution	NA

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