

Instruction Manual

Catalog Number	PK-CA707-99902
Description	<p>The human anticoagulant Annexin V is a 35-36 kDa, Ca²⁺-dependent phospholipid-binding protein with high affinity for phosphatidylserine (PS). During apoptosis, phosphatidylserine (PS) is translocated from the cytoplasmic face of the plasma membrane to the cell surface where it can be detected by labeled Annexin V which is a common probe for detecting apoptosis.</p> <p>5X Annexin V Binding Buffer is a concentrated stock solution used to prepare 1X Annexin V Binding Buffer which is optimized for staining cells with Annexin V conjugates.</p>
Quantity	15 ml
Excitation / Emission Maxima	NA
Molecular Structure	NA
Molecular Weight / Molecular Formula	NA
Purity	NA
Appearance / Formulation / Solubility	Colorless solution
Storage & Stability	Store at 4°C.
Applications	<p>Buffer optimized for binding assays with Annexin V conjugates, e.g. for use with the PromoKine Apoptotic/Necrotic Cells Detection Kit (Cat.# PK-CA707-30017) and Apoptotic/Necrotic/Healthy Cells Detection Kit (Cat.# PK-CA707-30018).</p> <p>PromoKine's 5X Annexin V Binding Buffer should be diluted to 1X with ultra-pure water before use. Rinse cells once with PBS, followed by 1X Annexin V Binding Buffer and incubate cells for 30 minutes in the dark. Rinse stained cells twice with 1X Annexin V Binding Buffer and visualize by fluorescence detection of choice.</p> <p>Staining concentrations for fluorescent Annexin V conjugates typically range from 0.25 – 2.5 µg/ml. We recommend determining the optimal staining concentration empirically. Generally, higher concentrations of Annexin V should be used for microscopy based assays and lower concentrations may be used for flow cytometry.</p> <p>For a complete Annexin V binding assay protocol, please refer to the protocols of the kits mentioned above.</p> <p>Note: Annexin V cannot be used to stain fixed cells or tissue sections.</p>
References	See kits PK-CA707-30017 and PK-CA707-30018.
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

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