

Instruction Manual

Catalog Number	PK-AB704-69006F
Quantity	100 tests
Clone	VAA-33
IgG Subtype	mouse IgG2ak
Purity	> 95% IgG2ak as determined by SDS-gel electrophoresis.
Specificity	Specifically recognizes Annexin V (VAC- α).
Cross-reactivity	There are no data about cross-reactivity with other species.
Formulation	50 mM TRIS 100 mM NaCl 1% BSA 0.02% NaN ₃ , pH 7.4
Storage	Stable at 2-8°C. After opening, aliquote contents and freeze at -20°C for long-term storage. Avoid repeated freeze/thaw cycles.
Shipping conditions	2-8°C
Applications	ELISA: PK-AB704-69006 is suitable for ELISA application. Western Blot: PK-AB704-69006 is suitable for Western blot analysis. Flow cytometry: PK-AB704-69006 is suitable as primary antibody for staining in FACS analysis. PK-AB704-69006F (FITC-labeled antibody) can be used for direct staining of cells. PK-AB704-69006B (biotinylated antibody) can be used in a two-step procedure using the biotin-(strept-)avidin system. Intracellular Staining: For intracellular stainings the use of PK-AB704-69006F is recommended. Optimal antibody concentrations should be determined by the individual laboratory for each application. This antibody is also available as biotinylated conjugate.

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