

Instruction Manual

Catalog Number	PK-AB704-69006
Quantity	100 µg
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	PBS
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	<p>ELISA: PK-AB704-69006 is suitable for ELISA application.</p> <p>Western Blot: PK-AB704-69006 is suitable for Western blot analysis.</p> <p>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. BPK-AB704-69006B (biotinylated antibody) can be used in a two-step procedure using the biotin-(strept-)avidin system.</p> <p>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.</p> <p>This antibody is also available as biotinylated or FITC-labeled conjugate.</p>

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