

Instruction Manual

Catalog Number	PK-AB704-69206
Quantity	100 µg
Clone	SM1/1
IgG Subtype	mouse IgG2a
Purity	> 95% IgG2a as determined by SDS-gel electrophoresis
Specificity	Specifically recognizes human APO-1.
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	Flow cytometry: PK-AB704-69206 is suitable as primary antibody in staining for FACS analysis. Functional studies: PK-AB704-69206 was found to induce apoptosis at 100 - 500 ng/ml in JURKAT-cells and SKW6.4 cells if secondary crosslinking with anti-mouse IgG is used. Induces apoptosis in human Fastransfectants without crosslinking. Optimal antibody concentrations should be determined by the individual laboratory for each application.

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