

Mouse Anti-Human Apoptosis Antigen 1/Fas (clone SM1/23, biotinylated)

Instruction Manual

Catalog Number	PK-AB704-69207B
Quantity	100 tests
Clone	SM1/23
IgG Subtype	mouse IgG2b
Purity	> 95% IgG2b as determined by SDS-gel electrophoresis.
Specificity	Specifically recognizes APO-1.
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	Functional studies: PK-AB704-69207 was found to block induction of apoptosis induced by clone SM1/1 or other commercial anti APO-1 antibodies. Flow cytometry: PK-AB704-69207 is suitable as primary antibody for staining in FACS analysis. PK-AB704-69207F can be used for direct staining of cells. PK-AB704-69207B (biotinylated antibody) can be used in a two-step procedure using the biotin-(strept-)avidin system. Optimal antibody concentrations should be determined by the individual laboratory for each application. This antibody is also available as FITC-labeled conjugate.

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