

# APO-1/Fas (SM1/23) mAb, FITC

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

## Instruction Manual

Catalog Number	PK-AB704-69207F
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 <sub>3</sub> , 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 (FITC-labeled antibody) 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 biotinylated conjugate.

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