

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

Instruction Manual

Catalog Number	PK-AB704-69208F
Quantity	100 tests
Clone	APO-1-1
IgG Subtype	mouse IgG1
Purity	> 95% IgG1 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	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	<p>Immunohistochemistry: APO-1-1 (PK-AB704-69208, IgG1) is an isotype-switch variant of the original clone APO-1-3 (IgG3). Since undesired side reactions in immunohistochemistry are minimal for IgG1 antibodies, PK-AB704-69208 is especially useful for immunohistology on cryostat sections and on cytospin preparations.</p> <p>Flow cytometry: PK-AB704-69208 is also suitable as primary antibody in staining for FACS analysis, PK-AB704-69208F (FITC-labeled antibody) can be used for direct staining of cells, PK-AB704-69208B (biotinylated antibody) can be used in a two-step procedure using the biotin-(strept-)avidin system. Recommended concentration for the unlabelled antibody is 1 - 2 µg/ml and a 1:20 dilution of the labelled antibodies, as determined on SKW6.4 cells.</p> <p>Optimal antibody concentrations should be determined by the individual laboratory for each application.</p> <p>This antibody is also available as biotinylated conjugate.</p>

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