

# MAP1 (IN) antibody (pAb)

## Rabbit Anti-Human/Mouse Modulator of Apoptosis 1

### Instruction Manual

Catalog Number	PK-AB718-3005
Synonyms	MAP1 Antibody:
Description	Apoptosis plays a major role in normal organism development, tissue homeostasis, and removal of damaged cells. Disruption of this process has been implicated in a variety of diseases such as cancer. Members of the Bcl-2 family are known to be critical regulators of this process. These proteins are characterized by the presence of several conserved motifs termed Bcl-2 homology (BH) domains. A related protein termed MAP-1 has recently been identified. This protein contains a BH3-like domain and induces caspase-dependent apoptosis in mammalian cells when overexpressed. It forms homodimers and associates with Bcl-2 family members such as Bax, Bcl-2, and Bcl-XL in vitro and in vivo. It has been suggested that MAP-1 associates with the tumor suppressor RASSF1A following death receptor activation, allowing a conformational change in Bax that leads to cellular apoptosis.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal MAP-1 antibody was raised against a 15 amino acid peptide from near the center of human MAP-1 (Genbank accession No. NP_000539).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. MAP1 antibody can be used for the detection of MAP-1 by Western blot at 1 to 4 µg/mL.
Images	Available upon request.
References	Lockshin RA, Osborne B, and Zakeri Z. Cell death in the third millennium. Cell Death Differ. 2000; 7:2-7. Cory S, Huang DCS, and Adams JM. The Bcl-2 family: roles in cell survival and oncogenesis. Oncogene 2003; 22:8590-607. Heiser D, Labi V, Erlacher M, et al. The Bcl-2 protein family and its role in the development of neoplastic disease. Exp. Geron. 2004; 39:1125-35. Tan KO, Tan KML, Chan S-L, et al. MAP-1, a novel proapoptotic protein containing a BH3-like motif that associates with Bax through its Bcl-2 homology domains. J. Biol. Chem. 2001; 276:2802-7.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3005P EL4 Cell Lysate, Cat. No. PK-AB718-1287 Bax Antibody (NT), Cat. No. PK-AB718-3351 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335

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