

Bmf (NT) antibody (pAb)

Rabbit Anti-Human/Mouse ATR (Anthrax toxin receptor, tumor endothelial marker 8, TEM8; IN)

Instruction Manual

Catalog Number	PK-AB718-3031
Synonyms	Bmf Antibody:
Description	Apoptosis is related to many diseases and development. Members in the Bcl-2 family are critical regulators of apoptosis by either inhibiting or promoting cell death. Bcl-2 homology 3 (BH3) domain is a potent death domain. BH3-only proteins, including Bad, Bid, Bik, Hrk, Bim, Noxa, and PUMA, form a growing subclass of the Bcl-2 family. A novel BH3-only protein was recently identified in human and mouse and designated Bmf (for Bcl-2-modifying factor). The BH3 domain in Bmf is required both for binding to Bcl-2 proteins and for triggering apoptosis. In healthy cells, Bmf associates with the dynein light chain 2 (DLC2) component of the myosin V motors and is sequestered by the cell's actin cytoskeleton. Disruption of the actin cytoskeleton, either by depolymerization of actin filaments or by detachment of cells from the extracellular matrix, triggers release and activation of Bmf, initiating the downstream apoptotic program. Bmf is constitutively expressed in many tissues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bmf antibody was raised against a synthetic peptide corresponding to 13 amino acids near the amino terminus of human Bmf (GenBank Accession number Q96LC9). This sequence differs from that of mouse by one amino acid.
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, ICC Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. Bmf antibody can be used for detection of Bmf by Western blot at 2 µg/mL. A band at approximately 25 kDa can be detected. Antibody can also be used for immunocytochemistry starting at 10 µg/mL.
Images	Available upon request.
References	Hunt A, Evan G. Apoptosis. Till death us do part. Science. 2001;293(5536):1784-5. Puthalakath H, Villunger A, O'Reilly LA, Beaumont JG, Coultas L, Cheney RE, Huang DC, Strasser A. Bmf: a proapoptotic BH3-only protein regulated by interaction with the myosin V actin motor complex, activated by anoikis. Science. 2001;293(5536):1829-32.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3031P; HepG2 Cell Lysate, Cat. No. PK-AB718-1204 Bmf Antibody (CT), Cat. No. PK-AB718-3029; Bim Antibody, Cat. No. PK-AB718-2065 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335; PUMA Antibody (CT), Cat. No. PK-AB718-3041

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