

Mcl-1 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Mcl-1 (Bcl-2 Protein Family Member)

Instruction Manual

Catalog Number	PK-AB718-3477
Synonyms	Mcl-1 Antibody: Myeloid cell leukemia 1
Description	Myeloid cell leukemia-1 (Mcl-1) is a member of the Bcl-2 family of proteins that can act to promote cell survival. While the mechanism by which Mcl-1 inhibits apoptosis is not known, it is thought that it may heterodimerize and neutralize pro-apoptotic members of the Bcl-2 family such as Bim or Bak. Mcl-1 was originally identified in differentiating myeloid cells, but has since been shown to be expressed in multiple cell types. Mcl-1 is essential for embryogenesis and for the development and maintenance of B and T lymphocytes in animals. Mcl-1 exists as at least two distinct isoforms designated Mcl-1L and Mcl-1S. In marked contrast to the larger isoform of Mcl-1, overexpression of Mcl-1S promotes cell death.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Mcl-1 antibody was raised against a peptide corresponding to 16 amino acids near the carboxy-terminus of human Mcl-1 (GenBank accession no. NP_068779).
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	Detects large isoform of Mcl-1 only
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, IF 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. Mcl-1 antibody can be used for detection of the large isoform of Mcl-1 only by Western blot at 1 to 2 µg/mL. Antibody can also be used for immunocytochemistry starting at 2 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Edwards SW, Derouet M, Howse M, et al. Regulation of neutrophil apoptosis by Mcl-1. <i>Biochem Soc Trans.</i> 2004; 32:489-92. Cuconati A, Mukherjee C, Perez D, et al. DNA damage response and MCL-1 destruction initiate apoptosis in adenovirus-infected cells. <i>Genes and Dev.</i> 2003; 17:2922-32. Kozopas KM, Yang T, Buchan HL, et al. MCL1, a gene expressed in programmed myeloid cell differentiation, has sequence similarity to BCL2. <i>Proc. Natl. Acad. Sci. USA</i> 1993; 90:3516-20.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3477P Raji Cell Lysate, Cat. No. PK-AB718-1207 Mcl-1 Antibody (NT), Cat. No. PK-AB718-3479; Mcl-1 Antibody (IN), Cat. No. PK-AB718-3481 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335 Bim Antibody (IN), Cat. No. PK-AB718-2065; Bak Antibody (NT), Cat. No. PK-AB718-3347

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