

OMI / HtrA2 (C2) antibody (pAb)

Rabbit Anti-Human OMI (IAP Inhibitor)

Instruction Manual

Catalog Number	PK-AB718-3051
Synonyms	OMI Antibody: HtrA2
Description	Inhibitor of apoptosis proteins (IAPs) were initially identified in baculoviruses as proteins that inhibit apoptosis of the host cells to allow time for viral replication. Cellular homologues containing at least one baculoviral IAP repeat (BIR) motif essential for their anti-apoptosis activity have been identified in yeasts and higher organisms and often act by binding and inhibiting processed caspases. The activity of these proteins can be modulated by the expression of proteins such as Smac/DIABLO and XAF-1 which displace or prevent the binding of caspases by IAPs. Recently, a mitochondrial serine protease termed Omi/HtrA2 has been found to bind IAPs. Similar to Smac, Omi possesses a conserved IAP-binding motif, but acts to cleave IAPs to irreversibly inactivate IAPs and promote apoptosis.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal OMI antibody was raised against a peptide corresponding to 16 amino acids near the C-terminus of human OMI (GenBank accession no. AAB94569).
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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, IHC, 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. OMI antibody can be used for detection of OMI by Western blot at 0.5 to 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Crook NE, Clem RJ, and Miller LK. An apoptosis inhibiting baculovirus gene with a zinc finger like motif. <i>J. Virol.</i> 1993; 67:2168-2174. Liston P, Fong WG, and Korneluk RG. The inhibitors of apoptosis: there is more to life than Bcl2. <i>Oncogene</i> 2003; 22:8568-80. Vaux DL and Silke J. Mammalian mitochondrial IAP binding proteins. <i>Biochem. Biophys. Res. Comm.</i> 2003; 304:499-504. Suzuki Y, Imai Y, Nakayama H, et al. A serine protease, HtrA2, is released from the mitochondria and interacts with XIAP, inducing cell death. <i>Mol. Cell</i> 2001; 8:613-21.
Images	Available upon request.
Related Products	OMI Blocking Peptide, Cat. No. PK-AB718-3051P U937 Lysate, Cat. No. PK-AB718-1215

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