

Instruction Manual

Catalog Number	PK-AB718-2479
Synonyms	MCG10 Antibody: MCG10, Poly rC binding protein 4, pcbp4, alpha-CP4
Description	MCG10 is a member of the KH-domain RNA-binding protein family. Proteins of this family also referred to as alpha-CPs, bind to RNA with specificity for C-rich pyrimidine regions. Alpha-CPs play important roles in post-transcriptional activities and have different cellular distributions. MCG10 is induced by p53 and DNA damage via two p53-responsive promoter elements. Its overexpression in transfected cells results in the suppression of cell proliferation by inducing apoptosis and cell cycle arrest in G2-M. Decreased expression of one isoform of MCG10 has been observed in many forms of lung cancer, leading to speculation that MCG10 can act as a tumor suppressor. Multiple alternatively spliced isoforms of MCG10 have been identified.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal MCG10 antibody was raised against a 17 amino acid peptide from near the carboxy terminus of human MCG10 (GenBank accession no. EAW65166).
Purification Method	Affinity chromatography purified via peptide column.
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, 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. MCG10 antibody can be used for detection of MCG10 by Western blot at 0.5 - 2 µg/mL. Antibody can also be used for immunocytochemistry starting at 1 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Makeyev AV and Liebhaber SA. The poly(C)-binding proteins: a multiplicity of functions and a search for mechanisms. RNA 2002; 8:265-78. Zhu J and Chen X. MCG10, a novel p53 target gene that encodes a KH domain RNA-binding protein, is capable of inducing apoptosis and cell cycle arrest in G2-M. Mol. Cell Biol. 2000; 20:5602-5618. Pio R, Zudaire I, Pino I, et al. a-CP-4, encoded by a putative tumor suppressor gene at 3p21, but not its alternative splice variant aCP-4a, is underexpressed in lung cancer. Cancer Res. 2004; 64:4171-9. (07-01D)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2479P HeLa Cell Lysate, Cat. No. PK-AB718-1201

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