

Instruction Manual

Catalog Number	PK-AB718-3677
Synonyms	ASK1 Antibody: Apoptosis signal-regulating kinase, MAPKKK, MEK kinase 5
Description	Mitogen-activated protein (MAP) kinase cascades are activated in response to various extracellular stimuli, including cytokines, growth factors and environmental stresses. A novel MAP kinase kinase (MAPKKK) was recently identified and designated ASK1 (for apoptosis signal-regulating kinase 1) and MAPKKK5. ASK1 activated two different subgroups of MAPKK, MKK4 and MKK6, which in turn activated c-Jun N-terminal kinase (JNK) and p38 MAP kinase, respectively. ASK1/MAPKKK5 is activated by TNFR and Fas through the interaction with members of the TRAF family and Fas-associated protein Daxx. Overexpression of ASK1 induced apoptotic cell death, and a catalytically inactive form of ASK1 inhibited TNF- α -induced apoptosis. ASK1 is expressed in variety of tissues and cell lines.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ASK1 antibody was raised against a 16 amino acid peptide from near the carboxy terminus of human ASK1 (Genbank accession No. BAA12684).
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 μ l 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. ASK1 antibody can be used for the detection of ASK1 by Western blot at 0.5 - 2 μ g/mL. Antibody can also be used for immunocytochemistry starting at 2 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Ichijo H, Nishida E, Irie K, et al. Induction of apoptosis by ASK1, a mammalian MAPKKK that activates SAPK/JNK and p38 signaling pathways. <i>Science</i> 1997; 275:90-4. Wang XS, Diener K, Jannuzzi D, et al. Molecular cloning and characterization of a novel protein kinase with a catalytic domain homologous to mitogen-activated protein kinase kinase kinase. <i>J. Biol. Chem.</i> 1996; 271:31607-11.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3677P SW1353 Cell Lysate, Cat. No. PK-AB718-1214 ASK1 Antibody (CT), Cat. No. PK-AB718-1151 ASK1 Antibody (NT), Cat. No. PK-AB718-3679
FOR IN VITRO RESEARCH ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	