

## Rabbit Anti-Human ASC (Novel CARD-Containing Adaptor Protein)

### Instruction Manual

Catalog Number	PK-AB718-2287
Synonyms	ASC Antibody:
Description	Apoptosis is regulated by death domain (DD) and/or caspase recruitment domain (CARD) containing molecules and a caspase family of proteases. CARD containing cell death regulators include RAIDD, RICK, BCL10, Apaf-1, ARC, caspase-9, and caspase-2. A novel CARD domain containing protein was recently identified in human and mouse and designated ASC and TMS1. Ectopic expression of ASC/TMS1 induced apoptosis through activation of caspase-9 and inhibited the survival of human breast cancer cells. Overexpression of ASC/TMS1 induced DNA fragmentation. ASC/TMS1 is expressed in a variety of human and mouse tissues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ASC antibody was raised against a synthetic peptide (RESQSYLVEDLERS) corresponding to amino acids 182 to 195 of human ASC .
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 <span style="float: right;">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.</span>
Images	ASC antibody can be used for detection of ASC/TMS1 by Western blot at 1 µg/mL. An approximately 25 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 5 µg/mL.
References	<p>Masumoto J, Taniguchi S, Ayukawa K, Sarvotham H, Kishino T, Niikawa N, Hidaka E, Katsuyama T, Higuchi T, Sagara J. ASC, a novel 22-kDa protein, aggregates during apoptosis of human promyelocytic leukemia HL-60 cells. <i>J Biol Chem.</i> 1999;274(48):33835-8.</p> <p>Masumoto J, Taniguchi S, Nakayama K, Ayukawa K, Sagara J. Murine Ortholog of ASC, a CARD-Containing Protein, Self-Associates and Exhibits Restricted Distribution in Developing Mouse Embryos. <i>Exp Cell Res.</i> 2001;262(2):128-133.</p> <p>Conway KE, McConnell BB, Bowring CE, Donald CD, Warren ST , Vertino PM. TMS1, a novel proapoptotic caspase recruitment domain protein, is a target of methylation-induced gene silencing in human breast cancers. <i>Cancer Res.</i> 2000;60(22):6236-42.</p> <p>McConnell BB, Vertino PM. Activation of a caspase-9-mediated apoptotic pathway by subcellular redistribution of the novel caspase recruitment domain protein TMS1. <i>Cancer Res.</i> 2000;60(22):6243-7. (WD0102)</p>
Images	Available upon request.
Related Products	ASC Blocking Peptide, Cat. No. PK-AB718-2287P HL60 Lysate, Cat. No. PK-AB718-1209

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