

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-3871
<b>Synonyms</b>	BAG-1 Antibody: Bcl-2-binding athanogene-1, receptor-associated protein 46, RAP46
<b>Description</b>	Bcl-2-associated athanogene 1 (BAG-1) was first identified as an anti-apoptotic bcl-2-binding protein. Later it was found to bind the molecular chaperones Hsp70 and Hsc70 through its carboxy-terminal sequence (termed the Bag domain), resulting in the inhibition of the refolding activity of these chaperones. It is thought that by binding and inhibiting these molecular chaperones, BAG-1 is able to modulate the expression level of proteins requiring chaperones to fold correctly. One such group of proteins that are affected is glucocorticoid receptors. Other reports have suggested that the level of BAG-1 expression correlates with the aggressiveness of various cancers. Multiple isoforms of BAG-1 are known to exist.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	BAG-1 antibody was raised against a 14 amino acid peptide from near the amino terminus of human BAG-1.
<b>Purification Method</b>	Affinity chromatography purified via peptide column.
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Human, Mouse
<b>Specificity</b>	
<b>Formulation</b>	Antibody is supplied in PBS containing 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	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.
<b>Applications</b>	E, WB, IHC, IF <b>Note:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.  BAG-1 antibody can be used for the detection of BAG-1 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2 µg/mL. For immunofluorescence start at 20 µg/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Takayama S, Sato T, Kraweski K, et al. Cloning and functional analysis of BAG-1: a novel Bcl2-binding protein with anti-cell death activity. <i>Cell</i> 1995; 80:279-84. Nollen EAA, Brunsting JF, Song J, et al. Bag1 functions in vivo as a negative regulator of Hsp70 chaperone activity. <i>Mol. Cell. Biol.</i> 2000; 20:1083-8. Cato ACB and Mink S. BAG-1 family of cochaperones in the modulation of nuclear receptor action. <i>J. Steroid Biochem. &amp; Mol. Biol.</i> 2001; 78:379-88. Kajewska M, Turner BC, Shabaik A, et al. Expression of BAG-1 protein correlates with aggressive behavior of prostate cancers. <i>Prostate</i> 2006; 66:801-10.
<b>Images</b>	Available upon request.
<b>Related Products</b>	BAG-1 Antibody, Cat. No. PK-AB718-3869 BAG-1 Peptide, Cat. No. PK-AB718-3871P Bcl-2Antibody, Cat. No. PK-AB718-3335 PC-3 Lysate, Cat. No. PK-AB718-1216

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