

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3869P
<b>Quantity</b>	50 µg
<b>Source</b>	14 amino acids near the carboxy terminus of human BAG1
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 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>	Store BAG-1 peptide at -20°C, stable for one year.
<b>Application</b>	BAG-1 peptide is used for blocking the activity of BAG1 antibody.
<b>References</b>	<p>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.</p> <p>Tsuneyoshi N, Fukudome K, Kohara J, et al. The functional and structural properties of MD-2 required for liposaccharide binding are absent in MD-1. <i>J. Immunol.</i> 2005; 174:340-4.</p> <p>Nagai Y, Kobayashi T, Motoi Y, et al. The radioprotective 105/MD-1 complex links TLR2 and TLR4/MD-2 in antibody response to microbial membranes. <i>J. Immunol.</i> 2005; 174:7043-9.</p>

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