

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3969P
<b>Quantity</b>	50 µg
<b>Source</b>	14 amino acids near the amino terminus of human sumo
<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 Sumo peptide at -20°C, stable for one year.
<b>Application</b>	Sumo peptide is used for blocking activity of sumo antibody.
<b>References</b>	Matunis MJ, Coutavas E, and Blobel G. A novel ubiquitin-like modification modulates the partitioning of the Ran-GTPase-activating protein RanGAP1 between the cytosol and the nuclear pore complex. <i>J. Cell Biol.</i> 1996; 135:1457-70.

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