

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4335P
<b>Quantity</b>	50 µg
<b>Source</b>	
<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 Strep-tag II peptide at -20°C, stable for one year.
<b>Application</b>	Strep-tag II peptide is used for blocking the activity of strep-tag II antibody.
<b>References</b>	Schmidt TGM, Koepe J, Frank R, et al. Molecular interactions between the Strep-tag affinity peptide and its cognate target, streptavidin. <i>J. Mol. Biol.</i> 1996; 255:753-6.

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