

Instruction Manual

Catalog Number	PK-AB718-4337P
Quantity	50 µg
Source	
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 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	Store Strep-tag II peptide at -20°C, stable for one year.
Application	Strep-tag II peptide is used for blocking the activity of strep-tag II antibody.
References	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.