

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-1125P
<b>Quantity</b>	50 µg
<b>Source</b>	17 amino acids near the carboxy terminus of human SIRP alpha
<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 SIRP alpha peptide at -20°C, stable for one year.
<b>Application</b>	SIRP alpha peptide is used for blocking the activity of SIRP alpha antibody.
<b>References</b>	<p>Kharitononkov A, et al. Nature 1997;386:181-186</p> <p>Marsters SA; Sheridan JP; Donahue CJ; Pitti RM; Gray CL; Goddard AD; Bauer KD; Ashkenazi A. Curr Biol, 1996;6:1669-76.</p> <p>Bodmer JL; Burns K; Schneider P; Hofmann K; Steiner V; Thome M; Bornand T; Hahne M; Schroter M; Becker K; et al. Immunity, 1997;6:79-88.</p>

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