

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4043P
<b>Quantity</b>	50 µg
<b>Source</b>	14 amino acids near the center of human SCO1
<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 SCO1 peptide at -20°C, stable for one year.
<b>Application</b>	SCO1 peptide is used for blocking the activity of SCO1 antibody.
<b>References</b>	Glerum DM, Shtanko A, and Tzagoloff A. SCO1 and SCO2 act as high copy suppressors of a mitochondrial copper recruitment defect in <i>Saccharomyces cerevisiae</i> . <i>J. Biol. Chem.</i> 1996; 271:20531-5.

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