

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4119P
<b>Quantity</b>	50 µg
<b>Source</b>	24 amino acids near the carboxy terminus of human STIM1
<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 STIM1 peptide at -20°C, stable for one year.
<b>Application</b>	STIM1 peptide is used for blocking the activity of STIM1 antibody.
<b>References</b>	Zhang SL, Yu Y, Roos J, et al. STIM1 is a Ca <sup>2+</sup> sensor that activates CRAC channels and migrates from the Ca <sup>2+</sup> store to the plasma membrane. Nature 2005; 437:902-5.

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