

Instruction Manual

Catalog Number	PK-AB718-3363P
Quantity	50 µg
Source	16 amino acids near the amino terminus of mouse ST2
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 ST2 peptide at -20°C, stable for one year.
Application	ST2 peptide is used for blocking the activity of the ST2 antibody.
References	<p>Lanahan A, Williams JB, Sanders LK, et al. Growth factor-induced delayed early response genes. <i>Mol Cell Biol.</i> 1992; 12:3919-29.</p> <p>Janeway CA Jr and Medzhitov R. Innate immune recognition. <i>Annu. Rev. Immunol.</i> 2002; 20:197-216.</p> <p>O'Neill LAJ, Fitzgerald FA, and Bowie AG. The Toll-IL-1 receptor adaptor family grows to five members. <i>Trends in Imm.</i> 2003; 24:286-9.</p> <p>Janeway CA Jr and Medzhitov R. Innate immune recognition. <i>Annu. Rev. Immunol.</i> 2002;20:197-216.</p>

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