

Instruction Manual

Catalog Number	PK-AB718-3367P
Quantity	50 µg
Source	15 amino acids near the carboxy terminus of mouse SIGIRR
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 SIGIRR peptide at -20°C, stable for one year.
Application	SIGIRR peptide is used for blocking the activity of the SIGIRR antibody.
References	<p>Thomassen E, Renshaw BR, and Sims JE. Identification and characterization of SIGIRR, a molecule representing a novel subtype of the IL-1R superfamily. <i>Cytokine</i> 1999; 11:389-99.</p> <p>Takeda K, Kaisho T, and Akira S. Toll-like receptors. <i>Annu. Rev. Immunol.</i> 2003;21:335-76.</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.