

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3841P
<b>Quantity</b>	50 µg
<b>Source</b>	14 amino acids near the carboxy terminus of human RP105
<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 RP105 peptide at -20°C, stable for one year.
<b>Application</b>	RP105 peptide is used for blocking the activity of RP105 antibody.
<b>References</b>	<p>Divanovic S, Trompette A, Atabani SF, et al. Inhibition of TLR-4/MD-2 signaling by RP105/MD-1. <i>J. Endotoxin Res.</i> 2005; 11:363-8.</p> <p>Meyers JH, Sabatos CA, Chakravarti S, et al. The TIM family regulates autoimmune and allergic diseases. <i>Trends Mol. Med.</i> 2005; 11:362-9.</p> <p>Meyers JH, Chakravarti S, Schlesinger D, et al. TIM-4 is the ligand for TIM-1, and the TIM-1-TIM4 interaction regulates T cell proliferation. <i>Nat. Immunol.</i> 2005; 6:455-64.</p>

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