

Instruction Manual

Catalog Number	PK-AB718-1123P
Quantity	50 µg
Source	15 amino acids near the carboxy terminus of human Eotaxin
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 Eotaxin peptide at -20°C, stable for one year.
Application	Eotaxin peptide is used for blocking the activity of eotaxin antibody.
References	<p>Kitaura M, Nakajima T, Imai T, et al. Molecular cloning of human eotaxin, an eosinophil-selective CC chemokine, and identification of a specific eosinophil eotaxin receptor, CC chemokine receptor 3. <i>J Biol Chem</i> 1996;271:7725-30.</p> <p>Hofmann K, Bucher P, Tschopp J. The CARD domain: a new apoptotic signalling motif. <i>Trends Biochem Sci</i> 1997;22:155-156 (RD1299)</p>

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