

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3587P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the carboxy terminus of human Precerebellin
<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 Precerebellin peptide at -20°C, stable for one year.
<b>Application</b>	Precerebellin peptide is used for blocking the activity of Precerebellin antibody.
<b>References</b>	<p>Urade Y, Oberdick J, Molinar-Rode R, et al. Precerebellin is a cerebellum-specific protein with similarity to the globular component C1q B chain. <i>Proc. Natl. Acad. Sci. USA</i> 1991; 88:1069-73.</p> <p>Navas-Nartin SR and Weiss S. Coronavirus replication and pathogenesis: Implications for the recent outbreak of severe acute respiratory syndrome (SARS), and the challenge for vaccine development. <i>J Neurovirol.</i> 2004;10:75-85.</p> <p>Arbely E, Khattari Z, Brotons G, et al. A highly unusual palindromic transmembrane helical hairpin formed by SARS coronavirus E protein. <i>J Mol. Biol.</i> 2004;3414:769-79.</p> <p>Shapiro L and Scherer PE. The crystal structure of a complement-1q family protein suggests an evolutionary link to tumor necrosis factor. <i>Curr. Biol.</i> 1998; 8:335-8.</p>

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