

Instruction Manual

Catalog Number	PK-AB718-3589P
Quantity	50 µg
Source	15 amino acids near the center of human Precerebellin
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 Precerebellin peptide at -20°C, stable for one year.
Application	Precerebellin peptide is used for blocking the activity of Precerebellin antibody.
References	Urade Y, Oberdick J, Molinar-Rode R, et al. Precerebellin is a cerebellum-specific protein with similarity to the globular component C1q B chain. Proc. Natl. Acad. Sci. USA 1991; 88:1069-73. Wong GW, Wang J, Hug C, et al. A family of Acrp30/ adiponectin structural and functional paralogs. Proc. Natl. Acad. Sci. USA 2004; 101:10302-7. Shapiro L and Scherer PE. The crystal structure of a complement-1q family protein suggests an evolutionary link to tumor necrosis factor. Curr. Biol. 1998; 8:335-8.

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