

Instruction Manual

Catalog Number	PK-AB718-3979P
Quantity	50 µg
Source	13 amino acids near the carboxy terminus of human PEN2
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 PEN2 peptide at -20°C, stable for one year.
Application	PEN2 peptide is used for blocking activity of PEN2 antibody.
References	Periz G and Fortini ME. Functional reconstitution of gamma-secretase through coordinated expression of presenilin, Nicastrin, aph-1, and pen-2. J. Neurosci. Res. 2004; 77:309-22.

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