

Instruction Manual

Catalog Number	PK-AB718-2279P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human F1A alpha
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 F1A alpha peptide at -20°C, stable for one year.
Application	F1A alpha peptide is used for blocking the activity of F1A alpha antibody.
References	<p>Chan SL, Tan KO, Zhang L, Yee KS, Ronca F, Chan MY, Yu VC. F1Aalpha, a death receptor-binding protein homologous to the Caenorhabditis elegans sex-determining protein, FEM-1, is a caspase substrate that mediates apoptosis. <i>J Biol Chem.</i> 1999;274(45):32461-8.</p> <p>Sinha S, et al. Purification and cloning of amyloid precursor protein beta-secretase from human brain. <i>Nature</i> 1999;402:537-40</p> <p>Yan R, et al. Membrane-anchored aspartyl protease with Alzheimer's disease beta-secretase activity. <i>Nature</i> 1999;402:533-7</p>

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