

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2247P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino amino terminus of human BACE2
<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 BACE2 peptide at -20°C, stable for one year.
<b>Application</b>	BACE2 peptide is used for blocking the activity of BACE2 antibody.
<b>References</b>	<p>Yan R, et al. Membrane-anchored aspartyl protease with Alzheimer's disease beta-secretase activity. <i>Nature</i> 1999;402:533-7.</p> <p>Kobayashi K, Hatano M, Otaki M, Ogasawara T, Tokuhisa T. Expression of a murine homologue of the inhibitor of apoptosis protein is related to cell proliferation. <i>Proc Natl Acad Sci USA</i> 1999;96:1457-62 (WD0600)</p>

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