

Instruction Manual

Catalog Number	PK-AB718-2283P
Quantity	50 µg
Source	14 amino acids of mouse RIP3
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 RIP3 peptide at -20°C, stable for one year.
Application	RIP3 peptide is used for blocking the activity of RIP3 antibody.
References	<p>Pazdernik NJ, Donner DB, Goebel MG, Harrington MA. Mouse receptor interacting protein 3 does not contain a caspase-recruiting or a death domain but induces apoptosis and activates NF-κB. <i>Mol Cell Biol.</i> 1999;19(10):6500-8.</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.