

Instruction Manual

Catalog Number	PK-AB718-2075P
Quantity	50 µg
Source	20 amino amino terminus of human RICK
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 RICK peptide at -20°C, stable for one year.
Application	RICK peptide is used for blocking the activity of RICK antibody.
References	<p>Inohara N, del Peso L, Koseki T, Chen S, Nunez G. RICK, a novel protein kinase containing a caspase recruitment domain, interacts with CLARP and regulates CD95-mediated apoptosis. <i>J Biol Chem</i> 1998;273:12296-300.</p> <p>Krieser RJ, Eastman A. The cloning and expression of human deoxyribonuclease II. A possible role in apoptosis. <i>J Biol Chem</i> 1998;273:30909-14</p> <p>Shiokawa D, Tanuma S. Cloning of cDNAs encoding porcine and human DNase II. <i>Biochem Biophys Res Commun</i> 1998;247:864-9</p>

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