

RICK / RIP2 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat RIP-Like Interacting CLARP Kinase

Instruction Manual

Catalog Number	PK-AB718-2075
Synonyms	RICK Antibody: RIP2
Description	Apoptosis is mediated by death domain (DD) and/or caspase recruitment domain (CARD) containing molecules and a caspase family of proteases. DD-containing serine/threonine kinase RIP regulates Fas-induced apoptosis. A novel CARD-containing serine/threonine kinase was recently identified and designated RICK/RIP2/CARDIAK for RIP-like interacting CLARP kinase, receptor interacting protein-2, and CARD-containing ICE associated kinase, respectively. RICK contains an N-terminal kinase catalytic domain and a C-terminal CARD domain. Overexpression of RICK induced apoptosis and activation of NF- κ B and JNK. RICK interacts with members of the TRAF family, CLARP and caspase-1. Thus, RICK represents a novel kinase that regulates TNF and Fas induced-apoptosis and that is involved in the generation of proinflammatory cytokine IL-1 β . The messenger RNA of RICK is expressed in multiple human tissues.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal RICK antibody was raised against a peptide corresponding to amino acids near the amino terminus of human RICK (Genbank accession no. O43353).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 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 μ l 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	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. RICK antibody can be used for detection of RICK by Western blot at 1 μ g/mL. An approximately 60 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 10 μ g/mL.
Images	Available upon request.
References	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. J Biol Chem 1998;273:12296-300 McCarthy JV, Ni J, Dixit VM. RIP2 is a novel NF- κ B-activating and cell death-inducing kinase. J Biol Chem 1998;273:16968-75 Thome M, Hofmann K, Burns K, Martinon F, Bodmer JL, Mattmann C, Tschopp J. Identification of CARDIAK, a RIP-like kinase that associates with caspase-1. Curr Biol 1998;8:885-8 (RD1299)
Images	Available upon request.
Related Products	A431 Lysate, Cat. No. PK-AB718-1202; RICK Peptide, Cat. No. PK-AB718-2075P Caspase-1 Antibody, Cat. No. PK-AB718-3459 RICK Antibody, Cat. No. PK-AB718-2183; RIP3 Antibody, Cat. No. PK-AB718-2283

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