

# RIP3 antibody (pAb)

## Rabbit Anti-Mouse/Rat Receptor Interacting Protein 3

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-2283
<b>Synonyms</b>	RIP3 Antibody:
<b>Description</b>	Certain serine/threonine protein kinases, such as ASK1, RIP, DAP, and ZIP kinases, are mediators of apoptosis. Receptor interacting proteins including RIP and RIP2/RICK mediate apoptosis induced by TNFR1 and Fas, two prototype members in the death receptor family. A novel member in the RIP kinase family was recently identified and designated RIP3. RIP3 contains N-terminal kinase domain but, unlike RIP or RIP2, lacks the C-terminal death or CARD domain. RIP3 binds to RIP and TNFR1, mediates TNFR1 induced apoptosis, and attenuates RIP and TNFR1 induced NF- $\kappa$ B activation. Overexpression of RIP3 induces apoptosis and NF- $\kappa$ B activation. The messenger RNA of RIP3 is expressed in a subset of adult tissues.
<b>Quantity</b>	100 $\mu$ g
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	Rabbit polyclonal RIP3 antibody was raised against a peptide (AQFGRGRGWQPFHK) corresponding to amino acids 473 to 486 of murine RIP3 .
<b>Purification Method</b>	Affinity chromatography purified via peptide column.
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Mouse, Rat
<b>Specificity</b>	
<b>Formulation</b>	Antibody is supplied in PBS containing 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 $\mu$ 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.
<b>Storage &amp; Stability</b>	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.
<b>Applications</b>	E, WB, IHC, IF 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. RIP3 antibody can be used for detection of RIP3 by Western blot at 1 $\mu$ g/mL. An approximately 57 kDa band can be detected. Antibody can also be used for immunohistochemistry starting at 5 $\mu$ g/mL. For immunofluorescence start at 20 $\mu$ g/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Yu PW, Huang BC, Shen M, Quast J, Chan E, Xu X, Nolan GP, Payan DG, Luo Y. Identification of RIP3, a RIP-like kinase that activates apoptosis and NF $\kappa$ B. <i>Curr Biol.</i> 1999;9(10):539-42. Sun X, Lee J, Navas T, Baldwin DT, Stewart TA, Dixit VM. RIP3, a novel apoptosis-inducing kinase. <i>J Biol Chem.</i> 1999;274(24):16871-5. 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- $\kappa$ B. <i>Mol Cell Bio.</i> 1999; 19(10):6500-8 (WD0102)
<b>Images</b>	Available upon request.
<b>Related Products</b>	3T3 (Balb) Lysate, Cat. No. PK-AB718-1212 RIP3 Peptide, Cat. No. PK-AB718-2283P

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