

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-1115
<b>Synonyms</b>	RAIDD Antibody: CRADD
<b>Description</b>	Apoptosis, or programmed cell death, occurs during normal cellular differentiation and development of multicellular organisms. Apoptosis is induced by certain cytokines including TNF and Fas ligand of the TNF family through their death domain (DD)-containing receptors, TNFR1 and Fas. The death signals are transduced by a group of DD-containing adapter molecules. A novel cell death adapter was recently identified by two independent groups and designated RAIDD (RIP-associated ICH-1/CED-3-homologous protein with DD) and CRADD (caspase and RIP adapter with DD). RAIDD contains a DD and a CARD (for caspase recruitment domain) which interact with RIP and caspase, respectively, to transduce death signals. RAIDD is constitutively expressed in many tissues and mediates apoptosis caused by Fas and TNFR-1.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	Rabbit polyclonal RAIDD antibody was raised against a peptide corresponding to amino acids 99 to 117 of human RAIDD.
<b>Purification Method</b>	Affinity chromatography purified via peptide column.
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Human
<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 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>	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, ICC, IF <span style="float: right;">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.</span> RAIDD antibody can be used for detection of RAIDD by Western blot at 0.5 - 1 µg/mL. A 22 kDa band should be detected. Antibody can also be used for immunocytochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Duan H, Dixit VM. RAIDD is a new 'death' adaptor molecule. <i>Nature</i> 1997;385:86-89 Ahmad M, Srinivasula SM, Wang L, Talanian RV, Litwack G, Fernandes-Alnemri T, Alnemri ES. CRADD, a novel human apoptotic adaptor molecule for caspase-2, and FasL/tumor necrosis factor receptor-interacting protein RIP. <i>Cancer Res</i> 1997 57:615-619 Hofmann K, Bucher P, Tschopp J. The CARD domain: a new apoptotic signalling motif. <i>Trends Biochem Sci</i> 1997;22:155-156 (RD1299)
<b>Images</b>	Available upon request.
<b>Related Products</b>	Caspase-2 Antibody, Cat. No. PK-AB718-3447 HeLa Lysate, Cat. No. PK-AB718-1201 RAIDD Antibody, Cat. No. PK-AB718-1117 RAIDD Peptide, Cat. No. PK-AB718-1115P

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