

Rabbit Anti-Human TNF Related Apoptosis Inducing Ligand (Ligand for DR4 & DR5)

Instruction Manual

Catalog Number	PK-AB718-1113
Synonyms	Trail Antibody: Apo-2L
Description	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 in the TNF family through their death domain containing receptors, TNFR1 and Fas. A novel member in the TNF family was recently identified and designated TRAIL (for TNF-related apoptosis-inducing ligand) and Apo-2L (for Apo-2 ligand). TRAIL is a type II membrane protein and expressed in a variety of human tissues. Two novel death domain containing receptors DR4 and DR5 have been identified as the receptor for TRAIL. Like TNF and Fas ligand, TRAIL induces apoptosis and NF- κ B activation in many tissues and cells.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal TRAIL antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human TRAIL .
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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, IHC !Note: 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. TRAIL antibody can be used for detection of TRAIL by Western blot at 1 μ g/mL dilution. Antibody can also be used for immunohistochemistry starting at 20 μ g/mL.
Images	Available upon request.
References	Wiley SR, Schooley K, Smolak PJ, et al. Identification and characterization of a new member of the TNF family that induces apoptosis. <i>Immunity</i> 1995; 3:673-682. Pitti RM, Marsters SA, Ruppert S, et al. Induction of apoptosis by Apo-2 ligand, a new member of the tumor necrosis factor cytokine family. <i>J. Biol. Chem.</i> 1996; 271:12687-90. Pan G, O'Rourke K, Chinnaiyan AM, et al. The receptor for the cytotoxic ligand TRAIL. <i>Science</i> 1997; 276:111-113. Schneider P, Thome M, Burns K, et al. TRAIL receptors 1 (DR4) and 2 (DR5) signal FADD-dependent apoptosis and activate NF- κ B. <i>Immunity</i> 1997; 7:831-836.
Images	Available upon request.
Related Products	DR4 Antibody (NT), Cat. No. PK-AB718-1166 HeLa Cell Lysate, Cat. No. PK-AB718-1201 DR4 Antibody (NT), Cat. No. PK-AB718-1166

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