

TL1A (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat TNF-like 1A (DR3 Ligand)

Instruction Manual

Catalog Number	PK-AB718-3109
Synonyms	TL1A Antibody; TL1A
Description	Members in the TNF and its receptor superfamilies regulate immune responses and induce apoptosis. DR3 (also termed Wsl-1, Apo-3, TRAMP, and LARD) is preferentially expressed by T lymphocytes and upregulated during T cell activation. The ligand for DR3 was recently identified and designated TL1A. TL1A also binds decoy receptor DcR3/TR6, which is expressed in several lung and colon carcinomas and in some normal tissues. TL1A induces apoptosis and NF- κ B activation in DR3 expressing cells, which is antagonized by DcR3. TL1A is upregulated by proinflammatory cytokines TNF and IL-1. TL1A is a longer variant of TL1 (also called VEGI).
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal TL1A antibody was raised against a peptide corresponding to 14 amino acids near the N-terminus of human TL1A (GenBank accession no. AAM77366).
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	This polyclonal antibody will not cross-react with TL1.
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 I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. TL1A antibody can be used for detection of TL1A by Western blot at 0.5 to 2 μ g/mL.
Images	Available upon request.
References	Chinnayan AM, O'Rourke K, Yu GL, et al. Signal transduction by DR3, a death domain-containing receptor related to TNFR-1 and CD95. <i>Science</i> 1996; 274:990-2. Kitson J, Raven T, Jiang YP, et al. A death-domain-containing receptor that mediates apoptosis. <i>Nature</i> 1996; 384:372-5. Screaton GR, Xu XN, Olsen AL, et al. LARD: a new lymphoid-specific death domain containing receptor regulated by pre-mRNA splicing. <i>Proc. Natl. Acad. Sci. USA</i> 1997; 94:4615-9. Bodmer JL, Burns K, Schneider P. TRAMP, a novel apoptosis-mediating receptor with sequence homology to tumor necrosis factor receptor 1 and Fas (Apo-1/CD95). <i>Immunity</i> 1997; 6:79-88.
Images	Available upon request.
Related Products	PC-3 Cell Lysate, Cat. No. PK-AB718-1216 TL1A Antibody, Cat. No. PK-AB718-3789. Dc3 Antibody, Cat. No. PK-AB718-2140. DR3 Antibody (NT), Cat. No. PK-AB718-1166.

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