

TCCR (NT) antibody (pAb)

Rabbit Anti-Human T-cell Cytokine Receptor

Instruction Manual

Catalog Number	PK-AB718-2481
Synonyms	TCCR Antibody: Type I T-cell cytokine receptor, IL-27 receptor alpha, IL-27Ra
Description	Upon antigen challenge, T-helper cells differentiate into two functional distinct subsets, Th1 and Th2. Th1 cells produce IL-2, IFN- γ and lymphotoxin- β that augment cell mediated immune response while Th2 cells secrete IL-4, IL-5, and IL-10 that enhance humoral immunity. The function of T-helper cells is regulated by cytokines. A novel cytokine receptor was recently identified and cloned. It is a new member in the type I cytokine receptor family and designated TCCR for T-cell cytokine receptor and WSX-1. TCCR deficient mice had impaired Th1 responses to protein antigen challenge, including decreased levels of IFN- γ and Th1-dependent antibody IgG2a. TCCR is predominately expressed in thymus, spleen, lymph nodes and peripheral blood leukocytes.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal TCCR antibody was raised against a synthetic peptide (GSAGPLQCYGVGPLGD) corresponding to amino acids 44 to 59 of human TCCR precursor .
Purification Method	Affinity chromatography purified via peptide column.
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, 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.
Images	TCCR antibody can be used for detection of TCCR by Western blot at 1 μ g/mL. An approximately 70 kDa band can be detected. Antibody can also be used for immunohistochemistry starting at 10 μ g/mL. For immunofluorescence start at 20 μ g/mL.
References	Chen Q, Ghilardi N, Wang H, Baker T, Xie MH, Gurney A, Grewal IS and de Sauvage FJ. Development of Th1-type immune responses requires the type I cytokine receptor TCCR Nature 2000;407(6806):916-920 Sprecher,C.A., Grant,F.J., Baumgartner,J.W., Presnell,S.R., Schrader,S.K., Yamagiwa,T., Whitmore,T.E., O'Hara,P.J. and Foster,D.F. Cloning and characterization of a novel class I cytokine receptor Biochem. Biophys. Res. Commun. 1998;246(1):82-90
Images	Available upon request.
Related Products	Spleen Lysate, Cat. No. PK-AB718-1306 TCCR Peptide, Cat. No. PK-AB718-2481P

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