

Carabin (NT) antibody (pAb)

Rabbit Anti-Human/Mouse Carabin (NT)

Instruction Manual

Catalog Number	PK-AB718-4269
Synonyms	Carabin Antibody; TBC1 domain family member 10C, TBC1D10C
Description	Antigen binding by the T-cell receptor (TCR) is one of the critical first steps in the immune response, triggering a cascade of signaling pathways that ultimately lead to T-cell activation. Screening a yeast two-hybrid screen of a human T-cell cDNA library with calcineurin, a protein phosphatase involved in multiple signaling pathways including T-cell activation, resulted in the identification of Carabin, a member of the TBC1 domain family of proteins, as a calcineurin-binding protein. Unlike other members of the TBC1 domain protein family which are thought to have a role in regulating cell growth and differentiation, further experiments demonstrated that Carabin is part of a negative regulatory loop for the intracellular TCR signaling pathway as well as an inhibitor of the Ras signaling pathway, suggesting that Carabin may also mediate crosstalk between calcineurin and Ras. Carabin antibody does not recognize TBC1D10A or TBC1D10B. Carabin is known to exist in multiple isoforms.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Carabin antibody was raised in rabbits against a 17 amino acid peptide from near the amino terminus of human Carabin.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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 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.
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. Carabin antibody can be used for detection of Carabin by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL.
Images	Available upon request.
References	Weil R and Israel A. Deciphering the pathway from the TCR to NF-kappaB. Cell Death Differ. 2006; 13:826-33. Im SH and Rao A. Activation and deactivation of gene expression Ca2+/calcineurin-NFAT-mediated signaling. Mol. Cells 2004; 18:1-9. Pan F, Sun L, Kardian DB, et al. Feedback inhibition of calcineurin and ras by a dual inhibitory protein carabin. Nature 2007; 445:433-6. White RA, Pasztor LM, Richardson PM, et al. The gene encoding TBC1D1 with homology to the tre-2/USP6 oncogene, Bub2, and cdc16 maps to mouse chromosome 5 and human chromosome 4. Cytogenet. Cell Genet. 2000; 89:272-5.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4269P; Carabin Peptide Cat.No. PK-AB718-1306; Human Spleen Tissue Lysate

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