

# CRTH2 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat CRTH2 (NT)

## Instruction Manual

Catalog Number	PK-AB718-4027
Synonyms	CRTH2 Antibody: Chemoattractant receptor-homologous molecule expressed on Th2 cells, GPR44
Description	The chemoattractant receptor-homologous molecule expressed on Th2 cells (CRTH2) is a recently identified receptor for the prostaglandin D2 (PGD2) in addition to the classic prostaglandin D receptor. CRTH2 is expressed on Th2 cells and eosinophils and mediates chemotaxis of these cells to PGD2 and is thus thought to be a key receptor mediating eosinophil and Th2 recruitment during allergic responses. However, CRTH2-null mice showed enhanced eosinophil recruitment into the lung consistent with observations that the CRTH2-null mice produced significantly higher amounts of interleukin-5 (IL-5) and IL-3. This suggests that CRTH2 plays a nonredundant role in restricting eosinophilia and allergic response in vivo. At least two different isoforms of CRTH2 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	CRTH2 antibody was raised in rabbits against a 18 amino acid peptide from near the amino terminus of human CRTH2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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, IF <b>INote:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. CRTH2 antibody can be used for detection of CRTH2 by Western blot a 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Nagata K, Hirai K, Tanaka K, et al. CRTH2, an orphan receptor of T-helper-2-cells, is expressed on basophils and eosinophils and responds to mast cell-derived factor(s). FEBS Lett. 1999; 459:195-9. Shichijo M, Sugimoto H, Nagao K, et al. Chemoattractant receptor-homologous molecule expressed on Th2 cells activation in vivo increases blood leukocyte counts and its blockage abrogates 13,14-dihydro-15-keto-prostaglandin D2-induced eosinophilia in rats. Chevalier E, Stock J, Fisher T, et al. Cutting edge: chemoattractant receptor-homologous molecule expressed on Th2 cells plays a restricting role on IL-5 production and eosinophil recruitment. J. Immunol. 2005; 2056-60. (07-01D)
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4027P; CRTH2 Peptide Cat.No. PK-AB718-1205; Jurkat Cell Lysate

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