

PRODUCT INFORMATION

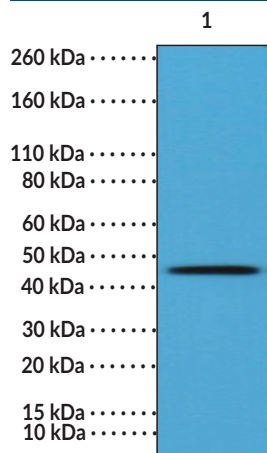


OX40/CD134 (C-Term) Rabbit Monoclonal Antibody (Clone RM313) Item No. 32255

Overview and Properties

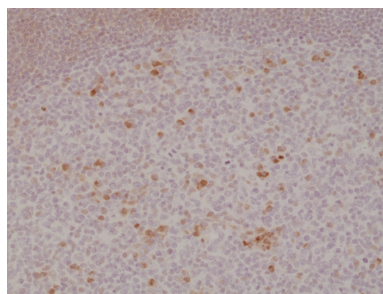
Contents:	This vial contains 100 µl protein A-affinity purified monoclonal antibody.
Immunogen:	Peptide from the C-terminal region of human OX40
Cross Reactivity:	(+) OX40
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	50% glycerol/PBS, with 1% BSA and 0.09% sodium azide
Clone:	RM313
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:200-1:500 for IHC and 1:1,000-1-2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Human spleen tissue lysate

WB of human spleen tissue lysate using OX40/CD134 (C-Term) Rabbit Monoclonal Antibody (Clone RM313) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using OX40/CD134 (C-Term) Rabbit Monoclonal Antibody (Clone RM313) at a 1:500 dilution.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/30/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

OX40, also known as TNF receptor superfamily member 4 (TNFRSF4), is a type I transmembrane glycoprotein and the cognate receptor for the T cell co-stimulatory molecule OX40L (TNFSF4; Item No. 32065).¹ OX40 assembles into trimers, with each monomer containing three cysteine-rich domains (CRDs) that bind OX40L and a cytoplasmic tail that interacts with TNF receptor-associated factor 1 (TRAF1), -2, -3, and -5 to induce a variety of intracellular signaling cascades, including NF- κ B, ERK, PI3K, and Akt pathways.^{1,2} OX40 is transiently expressed on CD4⁺ and CD8⁺ T cells in response to antigen stimulation in humans and is localized to lipid rafts.^{1,3} Binding of an OX40 trimer to trimeric OX40L, expressed by antigen-presenting cells (APCs), including dendritic cells (DCs), macrophages, and B cells, forms a hexameric OX40-OX40L complex that promotes effector T cell phenotype and function, proliferation, survival, and memory.³ Increased OX40 levels have been found in inflamed tissues isolated from patients with a variety of autoimmune diseases, including multiple sclerosis, systemic lupus erythematosus (SLE), or rheumatoid arthritis.³ Cayman's OX40 (C-Term) Rabbit Monoclonal Antibody (Clone RM313) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Buchan, S.L., Rogel, A., and Al-Shamkhani, A. The immunobiology of CD27 and OX40 and their potential as targets for cancer immunotherapy. *Blood* **131(1)**, 39-48 (2018).
2. Croft, M., So, T., Duan, W., *et al.* The significance of OX40 and OX40L to T-cell biology and immune disease. *Immunol. Rev.* **229(1)**, 173-191 (2009).
3. Webb, G.J., Hirschfield, G.M., and Lane, P.J.L. OX40, OX40L and autoimmunity: A comprehensive review. *Clin. Rev. Allergy Immunol.* **50(3)**, 312-332 (2016).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM