

# PRODUCT INFORMATION



## OX40/CD134 (human, recombinant)

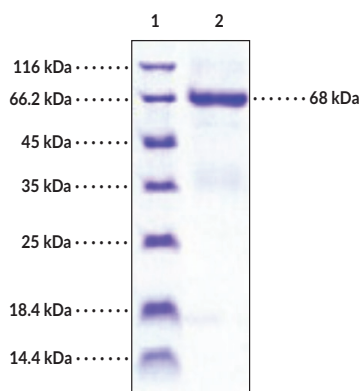
Item No. 31842

### Overview and Properties

<b>Synonyms:</b>	ACT35, IMD16, TNFRSF4, Tumor Necrosis Factor Receptor Superfamily Member 4 Precursor, TXGP1L
<b>Source:</b>	Active recombinant C-terminal human IgG1 Fc His-tagged OX40 expressed in HEK293 cells
<b>Amino Acids:</b>	29-216
<b>Uniprot No.:</b>	P43489
<b>Molecular Weight:</b>	48.2 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥85% estimated by SDS-PAGE
<b>Supplied in:</b>	Lyophilized from sterile PBS, pH 7.4
<b>Endotoxin Testing:</b>	<1.0 EU/μg, determined by the LAL endotoxin assay
<b>Protein Concentration:</b>	<i>batch specific</i> mg/ml
<b>Bioactivity:</b>	See figures for details

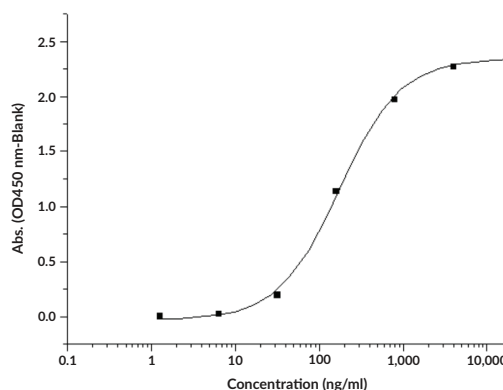
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: OX40/CD134

**SDS-PAGE Analysis of OX40/CD134.** This protein has a calculated molecular weight of 48.2 kDa. It has an apparent molecular weight of approximately 68 kDa by SDS-PAGE under reducing conditions due to glycosylation.



**OX40/CD134 Binding in Functional ELISA.** Immobilized cynomolgus mFc-TNFSF4 at 10 μg/ml (100 μl/well) can bind human OX40/CD134. The EC<sub>50</sub> value of human OX40/CD134 is 0.23-0.55 μg/ml.

**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, 04/22/2021

**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).<sup>1</sup> 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- $\kappa$ B, ERK, PI3K, and Akt pathways.<sup>1,2</sup> OX40 is transiently expressed on CD4<sup>+</sup> and CD8<sup>+</sup> T cells in response to antigen stimulation in humans and is localized to lipid rafts.<sup>1,3</sup> 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.<sup>3</sup> 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.<sup>3</sup> Cayman's OX40/CD134 (human, recombinant) protein can be used for binding assays. This protein is a disulfide-linked homodimer. The reduced monomer, composed of OX40 (amino acids 29-216) fused to His-tagged human IgG1 Fc at its C-terminus, consists of 436 amino acids, has a calculated molecular weight of 48.2 kDa, and a predicted N-terminus of Leu29 after signal peptide cleavage. As a result of glycosylation, the monomer migrates at approximately 68 kDa by SDS-PAGE under reducing conditions.

## 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