

CD134 (human):Fc (human), (recombinant)

The CD134 receptor (OX40) is a member of the tumor necrosis factor receptor (TNF-R) family. CD134 is a type I transmembrane receptor. It contains 3 extracellular cysteine-rich motifs. CD134 is mainly expressed in activated lymphocytes. After interaction with its ligand CD134L (OX40L), CD134 interacts with TNF receptor-associated factor 2 (TRAF2), TRAF5 and mediate activation of nuclear factors NF-κB and AP1. CD134-CD134L have been reported to be involved in T cell selection, migration and induction of Th1 and Th2 cytokine pattern. Moreover, it is involved in anti-tumor immunity, cell adhesion and various auto-immune diseases. Historical data has shown that CD134 receptor blocks T cell proliferation induced by CD134L.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-022-C050	50μg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

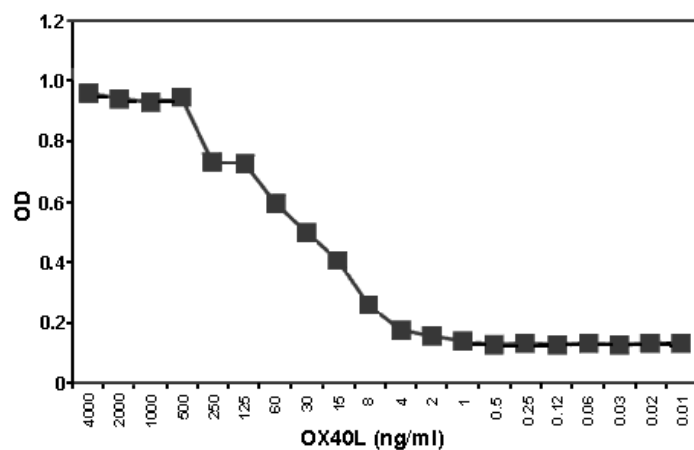


Figure: Receptor binding assay. Dose-dependent binding of recombinant human soluble CD134L to recombinant human CD134:Fc. Recombinant human CD134L saturates its receptor at a concentration range of 500-1'000ng/ml.**Method:** Receptor binding assay: A 96 well ELISA plate was coated over night with 50ng CD134 (human):Fc (human) (recombinant) (Prod. No. [ALX-522-022](#)) per well. After a blocking step, the indicated concentrations of CD134L, Soluble (human) (recombinant) (FLAG®) (Prod. No. ALX-522-023) were added for 1 hour. After extensive washing

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	OX40, TNFRSF 4
Application Notes	ELISA: binds to CD134L). Can be used together with anti-human IgG1 PAb to detect membrane-bound human CD134L. Has been used to ameliorate graft-versus-host disease, experimental allergic encephalomyelitis and inflammatory bowel disease.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~72kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The cysteine-rich region of human CD134 (OX40) (aa 29-214) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse CD134L (OX40L).
UniProt ID	P43489



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com