

CD137L (soluble) (human), (recombinant)

Biologically active CD137L (4-1BBL; TNFSF 9) comprising the whole extracellular tail domain (aa 49-93), which is important for the cross-linking of the protein that is essential for its biological activity.

The CD137 ligand (CD137L, 4-1BBL) is a member of the tumor necrosis factor (TNF) family. CD137L is a type II, transmembrane protein found on activated macrophages, dendritic cells and mature B cells. The interaction with its receptor 4-1BB induces recruitment of TNF receptor-associated factor 1 (TRAF1) and TRAF2 and interaction with the kinase p56^{lck}. CD137 and CD137L have been reported to be involved in tumor rejection, apoptosis, anti-viral immunity, diabetes, in T and B cell co-stimulation and modulation of the immune response.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-111-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

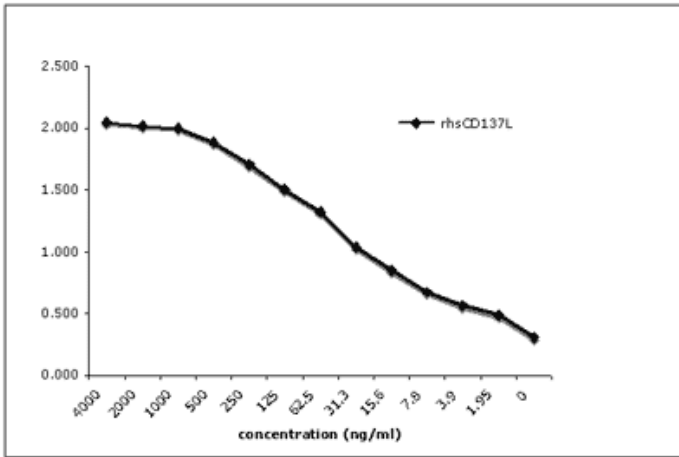


Figure: Dose-dependent binding of recombinant human soluble CD137L to CD137:Fc (human). Recombinant human CD137L saturates its receptor at a concentration range of 2-4'000ng/ml. **Method:** Receptor binding assay: a 96-well ELISA plate was coated O/N with 50ng per well of CD137 (human):Fc (human) (rec.) (4-1BB:Fc) (Prod. No. [ALX-522-031](#)). After a blocking step, the indicated concentrations of rhsCD137L were added for 1 hour. After extensive washing, Enhancer for Ligands (Prod. No. ALX-804-034) was added for 1 hour. Bound ligand was revealed by a rabbit anti-mouse IgG-HRP (1:1'000 dilution)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, 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	4-1BBL, TNFSF 9
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	≤1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS, 5% trehalose, and 5% mannitol.
MW	~27kDa
Purity	≥85% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl sterile distilled water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in HEK 293 cells. The extracellular domain of human CD137L (aa 49-254) is fused at the N-terminus to a linker peptide (10 aa) and a FLAG [®] -tag.
Specificity	Binds to human CD137 (4-1BB).
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

P41274



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

Last modified: May 29, 2024