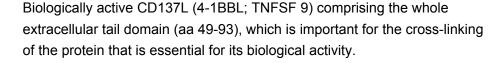
CD137L (soluble) (human), (recombinant)



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
ALX-522-111-3010	SuperPack - 3x10µg

Manuals, SDS & CofA

View Online »

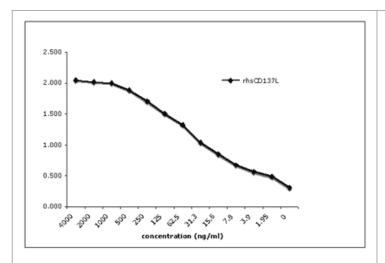


Figure: Dose-dependent binding of recombinant human soluble CD137L to CD137:Fc (human). Recombinant human CD137L saturates its receptor at a concentration range of 2-4'000ng/ml. **Method:** Receptor binding assay: a 96-well ELISA plate was coated O/N with 50ng per well of CD137 (human):Fc (human) (rec.) (4-1BB:Fc) (Prod. No. <u>ALX-522-031</u>). 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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

P41274

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

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

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