

CD40L, soluble (mouse) (recombinant) set

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-075-KI01	1Set
------------------	------

Manuals, SDS & CofA

[View Online »](#)

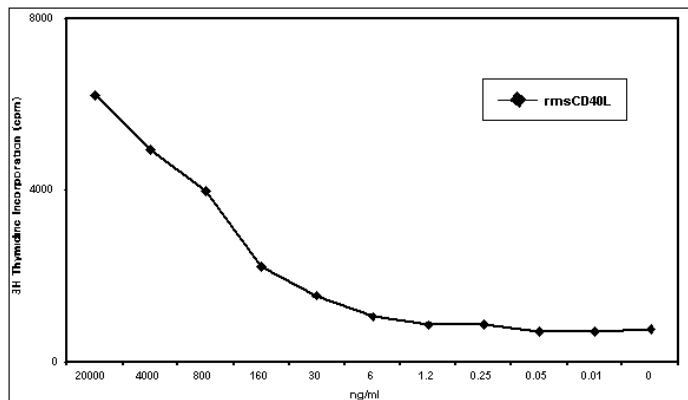


Figure: CD40L-induced proliferation of murine B cells.

Method: Splenocytes from Balb/c mice were purified on B220 magnetic beads using the MACS system. The resulting B cells were put in culture at a density of 70'000 cells/well in a 96-well plate. Cells were activated with the indicated concentrations of CD40L, Soluble (mouse) (recombinant) (Prod. No. [ALX-522-070](#)) in the presence of 1µg/ml of Enhancer for Ligands (Prod. No. ALX-804-034). After 36 hours

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	CD154, TNFSF 5, gp39
Application	ELISA, FUNC
Application Notes	Functional Application: Stimulates the proliferation of mouse B cells and dendritic cells. The activity of CD40L increases 1'000-fold (stimulation in the ng/ml range) in the presence of the cross-linking enhancer (Prod. No. ALX-804-034).
Concentration	CD40L: 0.1mg/ml after reconstitution.Enhancer: 1mg/ml after reconstitution.
Contents	10µg of CD40L, Soluble (mouse) (recombinant) (Prod. No. ALX-522-070) and 2×50µg of Enhancer for Ligands (Prod. No. ALX-804-034) which increases the biological activity of CD40L at least 1'000-fold.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
Purity	≥90% (SDS-PAGE).
Reconstitution	Reconstitute CD40L with 100µl sterile water and each vial of enhancer with 50µl of sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of mouse CD40L (CD154) (aa 115-260) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Specificity	CD40L binds to human and mouse CD40 in an ELISA assay.

Technical Info / Product Notes

CD40L (Prod. No. ALX-522-070) and Enhancer for Ligands (Prod. No. ALX-804-034) should be incubated together for 30 min at room temperature prior to cell application. Effects are typically best seen with 0.5-10µg/ml of CD40L together with 1-2µg/ml of enhancer. This may vary by cell type and experimental conditions. FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

P27548 (CD40L)



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)