

CD40L, soluble (human) (recombinant) set

Ligand plus enhancer for improved stability
and enhanced immune activation.

Citations: 19

[View Online »](#)

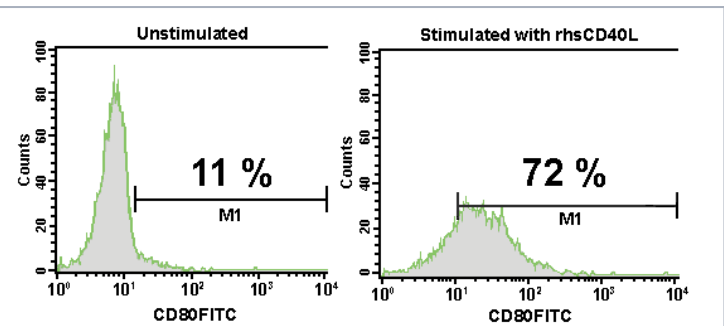
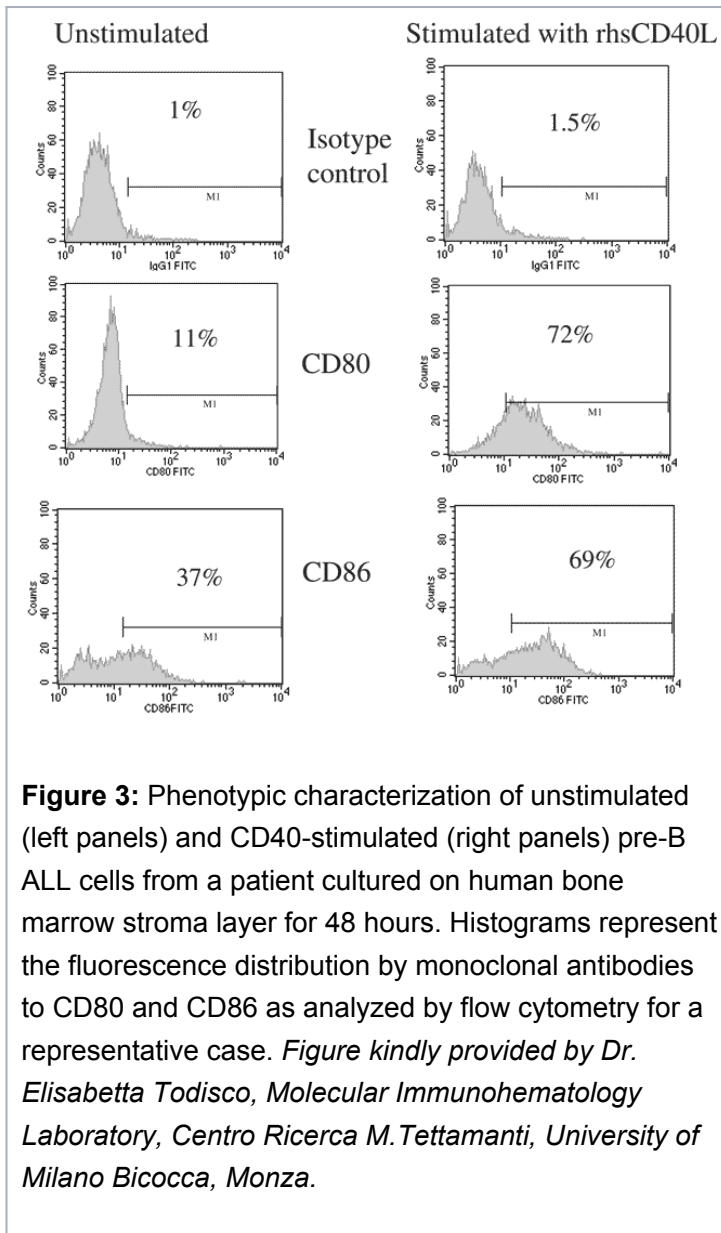
Ordering Information

[Order Online »](#)

ALX-850-064-KI01	1Set
------------------	------

Manuals, SDS & CofA

[View Online »](#)



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
Application	FUNC
Application Notes	Can be used to activate human B cells and dendritic cells. Stimulates growth of human B cells. For stimulation of mouse cells via CD40 use MAb to CD40 (mouse) (FGK45) (Prod. No. ALX-805-046).
Concentration	CD40L: 0.1mg/ml after reconstitutionEnhancer: 1mg/ml
Contents	10µg of CD40L, Soluble (human) (recombinant) (Prod. No. ALX-522-015) and 2×50µg of Enhancer for Ligands (Prod. No. ALX-804-034), which increases the biological activity of rhsCD40L at least 1'000-fold.
Formulation	CD40L: Lyophilized. Contains PBS. Enhancer: Liquid. In PBS, pH 7.4, containing 50% glycerol. The formulation contains no antimicrobial preservatives.
Purity	≥95% (SDS-PAGE).
Reconstitution	Reconstitute CD40L with 100µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or carrier protein.
Source	Produced in <i>E. coli</i> . The extracellular domain of human CD40L (CD154) (aa 116-261) is fused at the N-terminus to a linker peptide (6 aa) and a FLAG [®] -tag.
Specificity	CD40L binds to human CD40.

**Technical Info / Product
Notes**

CD40L (Prod. No. ALX-522-015) 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

P29965



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)