

Fc (human):CD40L (soluble) (human), (recombinant)

Ordering Information

[Order Online »](#)

ALX-522-076-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

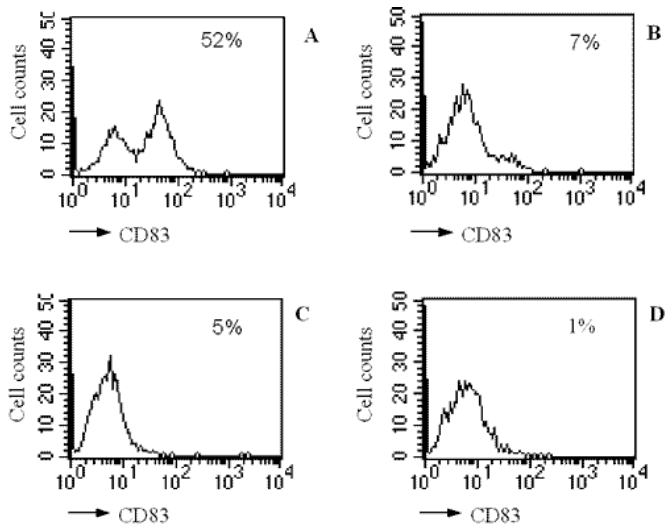


Figure: Proliferation of human dendritic cells (DC) induced by Fc (human):CD40L, Soluble (human) (rec.) (Prod. No. [ALX-522-076](#)). **Method:** Human DC were derived from GM-CSF/IL-4 treated monocytes. The so obtained DC, 10^5 cells/assay, were incubated with 3 μ g/ml Fc (human):CD40L, Soluble (human) (rec.) (A), 3 μ g/ml boiled Fc (human):CD40L, Soluble (human) (rec.) (B), 3 μ g/ml control:Fc (C) and media (D). After 18 hours, DC maturation was analyzed by immunostaining with a CD83-specific MAb and cytofluorimetry analysis. Indicated are the percentages of activated DC. "''", Fc (human):CD40L (soluble) (human)

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 Notes	ELISA: binds to human CD40:COMP at 2-50 ng/ml.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; BioWhittaker).
Formulation	Lyophilized. Contains PBS.
MW	~58kDa (SDS-PAGE).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl 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 extracellular domain of human CD40L (CD154) (aa 116-261) is fused at the N-terminus to the Fc portion of human IgG1 and a linker peptide (20 aa).
Specificity	Binds to human CD40.
Technical Info / Product Notes	Historical data has shown that Fc (human):CD40L stimulates the proliferation of human dendritic cells. Note: Does not require an enhancer.



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

P29965
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