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

Ordering Information

Order Online »

ALX-522-076-C010

10µg

Manuals, SDS & CofA

View Online »

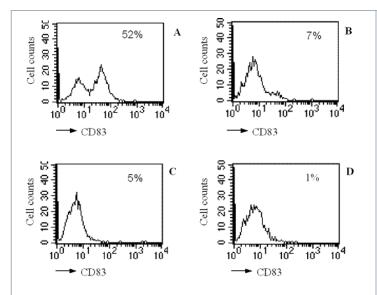


Figure: Proliferation of human dendritic cells (DC) induced by Fc (human):CD40L, Soluble (human) (rec.) (Prod. No. ALX-522-076).**Method:** Human DC were derived from GMCSF/IL-4 treated monocytes. The so obtained DC, 10⁵ cells/assay, were incubated with 3μg/ml Fc (human):CD40L, Soluble (human) (rec.) (A), 3μg/ml boiled Fc (human):CD40L, Soluble (human) (rec.) (B), 3μg/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.

UniProt ID P29965

Last modified: May 29, 2024



info-

eu@enzolifesciences.com