

# LAG-3 (human):Fc (human), (recombinant)

The lymphocyte activation gene-3 (LAG-3, CD223), a member of the immunoglobulin superfamily (IgSF) related to CD4, binds to the major histocompatibility complex (MHC) class II molecules but with higher affinity than CD4. Several alternative mRNA splice-variants of human LAG-3 have been described, two of them encoding potential secreted forms: LAG-3V1 (i.e. the D1-D2 domains of the protein, 36 kDa) and LAG-3V3 (D1-D3, 52 kDa). The longer form was detected by ELISA in the serum of healthy individuals as well as of tuberculosis patients with a favorable outcome. LAG-3 expression by T cell clones correlated with IFN- $\gamma$  production, and hence soluble LAG-3 has been suggested as a serological marker of Th1 responses.

Citations: 7

[View Online »](#)

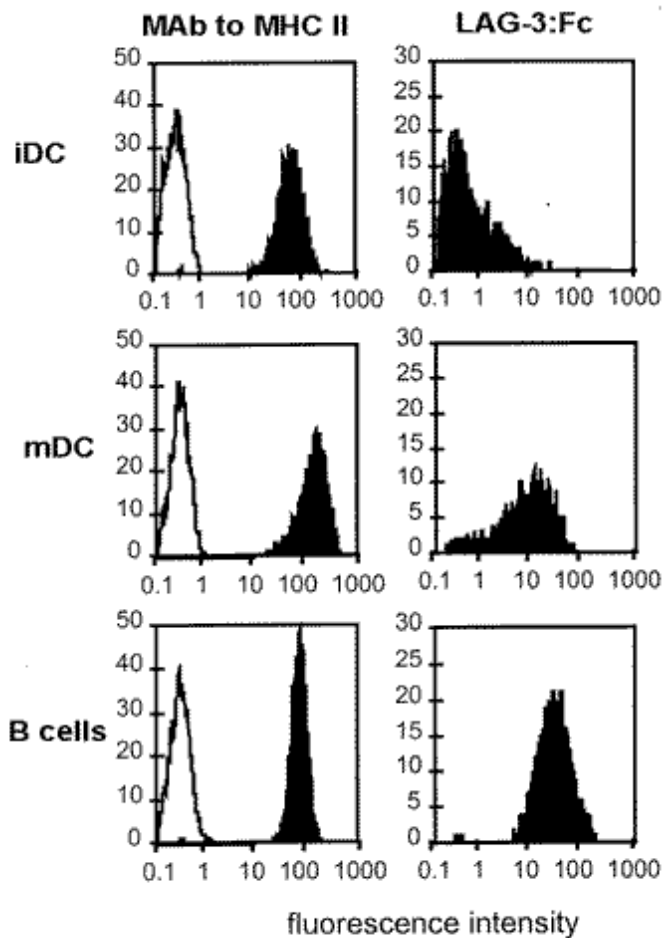
## Ordering Information

[Order Online »](#)

ALX-522-078-C050	50 $\mu$ g
------------------	------------

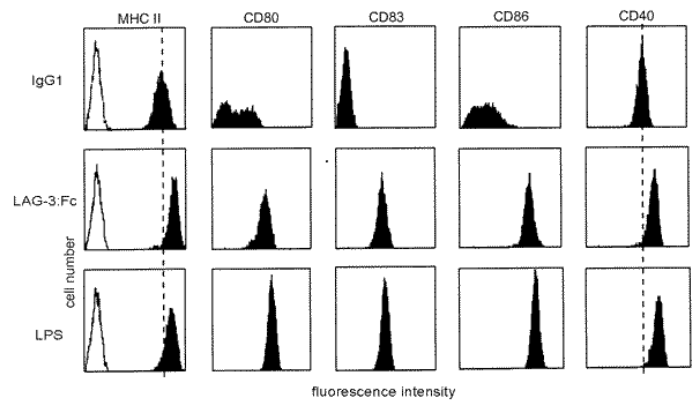
## Manuals, SDS & CofA

[View Online »](#)



**Figure 1:** Binding of LAG-3 (human):Fc (human) (rec.) (Prod. No. [ALX-522-078](#)) on dendritic and B cells.

**Method:** Monocyte-derived immature DC (iDC), DC matured for 48h with LPS (mDC) and EBV-transformed B-cells are incubated with either an anti-MHC class II MAb (10µg/ml) or LAG-3 (human):Fc (human) (rec.) (10µg/ml) for 30 min. at 4°C in PBS/1% BSA. The cells are then stained for 30 min. at 4°C with FITC-labeled goat anti-human Fc antibodies. The open histograms correspond negative control labeling [3]. " ", LAG-3 (human):Fc (human)



**Figure 2:** rhLAG-3:Fc-induced dendritic cells (DC) maturation. **Method:** Immature DC derived from human monocytes, resuspended at  $1 \times 10^6$  cells/ml in complete culture medium, are incubated with LAG-3 (human):Fc (human) (rec.) (Prod. No. [ALX-522-078](#)) (10µg/ml), LPS (positive control, 5µg/ml) or human IgG1 (isotype-matched negative control, 10µg/ml) for 48 hours. DC maturation induced by rhLAG-3:Fc binding to MHC class II or LPS is assessed by the increased expression of MHC Class II, CD80, CD83, CD86 and CD40 [3]. " ", LAG-3 (human):Fc (human)

## Handling & Storage

**Use/Stability** Stable for at least 6 months after receipt when stored at -20°C.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Lymphocyte activation gene-3, FDC protein, CD223

**Application Notes** Flow Cytometry: Detection of MHC class II molecules in combination with fluorescently labelled Ab to IgG1.

**Biological Activity** Inhibits binding of MAb to LAG-3 (human) to LAG-3. Induces maturation of human dendritic cells.

**Endotoxin Content** <0.1EU/mg purified protein (LAL test; Bio Whittaker).

**Formulation** Liquid. In PBS.

**MW** ~80kDa (SDS-PAGE).

**Purity** ≥99% (SDS-PAGE)

**Source** Produced in CHO cells. The sequence coding for the 4 extracellular Ig-like domains of human LAG-3 (D1-D4) is fused to the Fc portion of human IgG1.

**Specificity** Binds to human, mouse and monkey MHC class II.

**UniProt ID** P18627



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)