## 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-γ production, and hence soluble LAG-3 has been suggested as a serological marker of Th1 responses.

Citations: 7

**View Online »** 

**Ordering Information** 

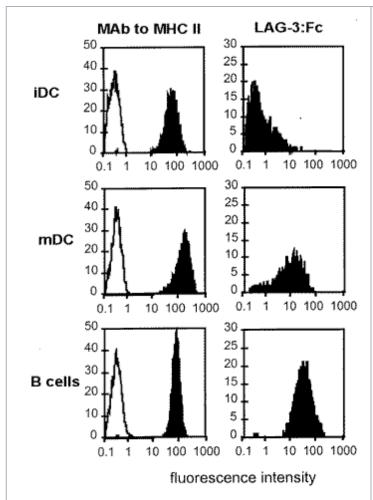
**Order Online** »

ALX-522-078-C050

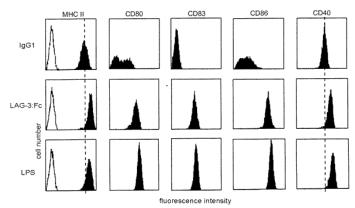
50µg

Manuals, SDS & CofA

**View Online** »



**Figure 1:** Binding of LAG-3 (human):Fc (human) (rec.) (Prod. No. <u>ALX-522-078</u>) 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 1×10<sup>6</sup> cells/ml in complete culture medium, are incubated with LAG-3 (human):Fc (human) (rec.) (Prod. No. <u>ALX-522-078</u>) (10μg/ml), LPS (positive control, 5μg/ml) or human lgG1 (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

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com