

# LAG-3 (mouse):Fc (mouse), (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: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-522-099-C050	50 $\mu$ g
------------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## 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

**Biological Activity** Induces maturation of mouse dendritic cells *in vitro*.  
Induces CD4 and CD8 antigen-specific T cells responses when coinjected s.c. in mice, together with the antigen.

**Concentration** 0.5mg/ml

**Endotoxin Content** <1EU/μg purified protein (LAL test; Bio Whittaker).

**Formulation** Liquid. In PBS.

**MW** ~180kDa (SPS-PAGE, nonreducing), ~80kDa (SPS-PAGE, reducing)

**Purity** ≥95% (SDS-PAGE)

**Reconstitution** Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.

**Source** Produced in CHO cells. The sequence coding for the 4 extracellular Ig-like domains of mouse LAG-3 (D1-D4) is fused to the Fc portion of mouse IgG2a (hinge-CH<sub>2</sub>-CH<sub>3</sub>).

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

**UniProt ID** Q61790



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 949 946 9120  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)