

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- γ 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 μ 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-CH₂-CH₃).

Specificity Binds to human and mouse MHC class II.

UniProt ID Q61790



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)