

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 »](#)

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 <i>in vitro</i> . 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 ₂ -CH ₃).
Specificity	Binds to human and mouse MHC class II.
UniProt ID	Q61790



ENZO LIFE SCIENCES,
INC.
Phone: +1 949 948 9120
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

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

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

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