

LAG-3 (human) recombinant monoclonal antibody (L4-PL33) (FITC conjugate)

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.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

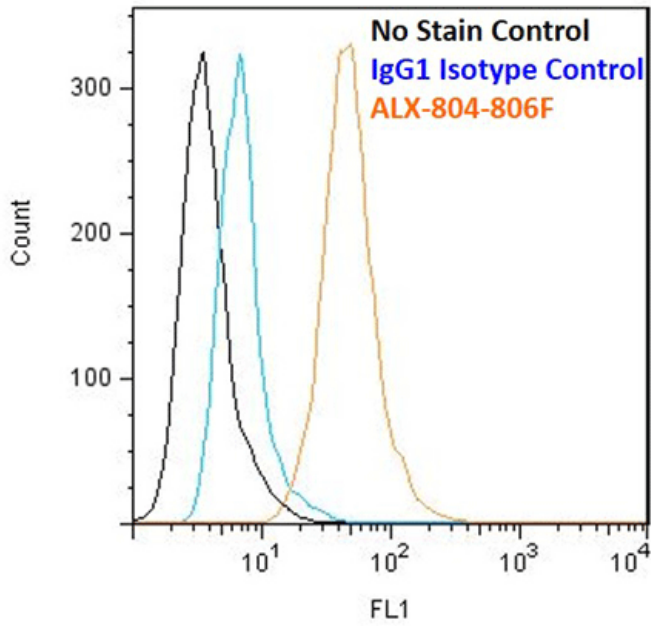
[Order Online »](#)

ENZ-ABS677F-0100

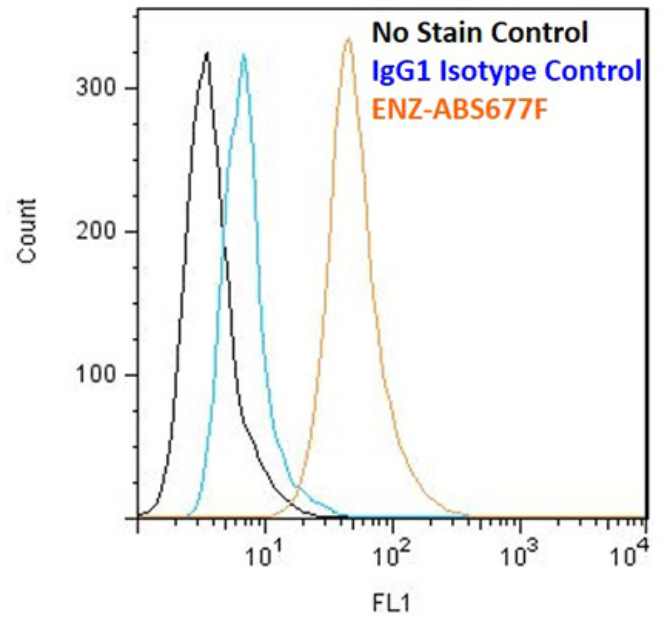
100 μ g

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 106 HDLM-2 cells stained using LAG-3 mAb (L4-PL33) (FITC conjugate), (Prod. No. ENZ-ABS677F), at a concentration of 10µg/ml, side-by-side comparison to LAG-3 mAb (17B4) (FITC conjugate), (Prod. No. ALX-804-806F).



Flow cytometry analysis of 106 HDLM-2 cells stained using LAG-3 mAb (L4-PL33) (FITC conjugate), (Prod. No. ENZ-ABS677F), at a concentration of 10µg/ml, side-by-side comparison to LAG-3 mAb (17B4) (FITC conjugate), (Prod. No. ALX-804-806F).

Handling & Storage

Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Lymphocyte activation gene-3, FDC protein, CD223
Application	Flow Cytometry
Clone	L4-PL33
Formulation	Liquid. In PBS containing 50mM Tris-HCl, 1% BSA and 0.02% sodium azide.
GenBank ID	NM_002286
Host	Mouse
Immunogen	Synthetic peptide corresponding to 30 aa (GPPAAAPGHPLAPGPHPAAPSSWGPRPRRY) from the first N-terminal D1 domain of human LAG-3 (lymphocyte activation gene-3).
Isotype	IgG1k
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from tissue culture supernatant.
Species Reactivity	Human

Specificity

Recognizes the 30 aa extra-loop of the first N-terminal D1 domain of LAG-3.

UniProt ID

P18627

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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)