LAG-3 (human) recombinant monoclonal antibody (L4-PL33) (PBS only)

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

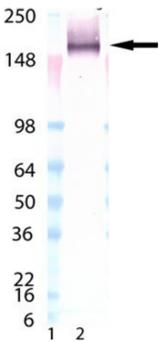
Order Online »

ENZ-ABS677PB-0100

100µg

Manuals, SDS & CofA

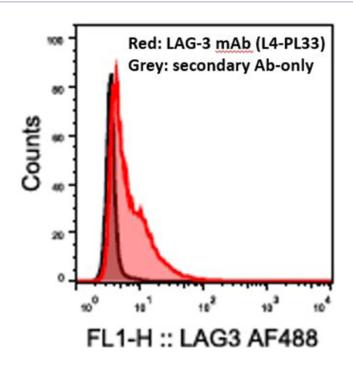
View Online »





Western blot analysis of LAG-3 mAb (L4-PL33) (Prod. No. ENZ-ABS677): Lane 1: MW marker, Lane 2: LAG-3 (human):Fc (human), (recombinant) (Prod. No. ALX-522-078).

Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human tonsil tissue stained with LAG-3 (human) recombinant mAb (L4-PL33) at 6µg/ml.



Flow cytometry analysis of 0.5×10⁶ THP-1 cells stained using LAG-3 mAb (L4-PL33) at a concentration of 5µg/ml, followed by Alexa Fluor® 488 conjugated anti-mouse secondary antibody.

Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored as recommended.

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°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 ELISA, Flow Cytometry, ICC, IHC, IP, WB

Clone L4-PL33

Formulation Liquid. In PBS-only. Preservative and stabilizer free.

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 IgG1ĸ

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:100)Immunohistochemistry (frozen

sections, 1:150; FFPE sections, 1:150)Western Blot (1-5

 μ g/ml)Immunoprecipitation (10 μ g/ml)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,

Phone: 800.942.0430

info-