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

Geomptote Stizt b 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

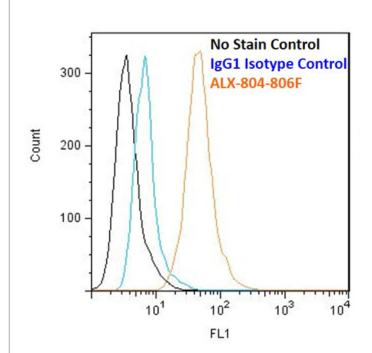
Order Online »

ENZ-ABS677F-0100

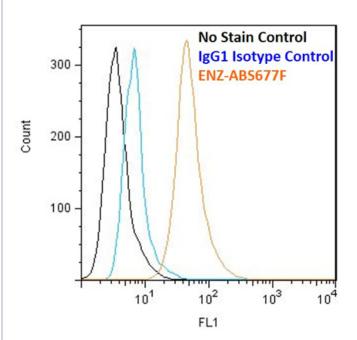
100µg

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 NameLymphocyte 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 IgG1ĸ

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



info-