## LAG-3 (human) recombinant monoclonal antibody (L4-PL33) (ATTO647N

General depends (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** 

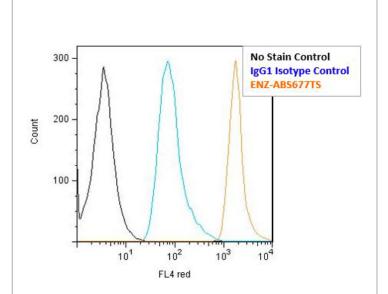
Order Online »

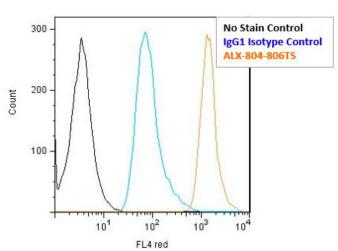
ENZ-ABS677TS-0100

100µg

Manuals, SDS & CofA

**View Online** »





Flow cytometry analysis of 10<sup>6</sup> HDLM-2 cells stained using LAG-3 mAb (L4-PL33) (ATTO 647N conjugate), (Prod. No. ENZ-ABS677TS), at a concentration of 10μg/ml, side-by-side comparison to LAG-3 mAb (17B4) (ATTO 647N conjugate), (Prod. No. ALX-804-806TS).

Flow cytometry analysis of  $10^6$  HDLM-2 cells stained using LAG-3 mAb (L4-PL33) (ATTO 647N conjugate), (Prod. No. ENZ-ABS677TS), at a concentration of  $10\mu g/ml$ , side-by-side comparison to LAG-3 mAb (17B4) (ATTO 647N conjugate), (Prod. No. ALX-804-806TS).

## **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, ICC

Clone 17B4

Formulation Liquid. In PBS containing 50mM Tris, 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).

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 hybridoma 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-