LAG-3 (human) monoclonal antibody (17B4) (preservative free)

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.

Citations: 9

View Online »

Ordering Information

Order Online »

ALX-804-806PF-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored at -20°C.

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 Flow Cytometry, FUNC, ICC, IHC, IP, WB

Application Notes Functional Application: Inhibition

Clone 17B4

Formulation Liquid. In PBS containing 10% glycerol.

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).

lgG1

Purity Detail Purified from concentrated hybridoma tissue culture

supernatant by Protein A Sepharose™ CL-4B Flow Fast

Column.

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections,

1:150)Immunoprecipitation (10µg/ml)Western Blot (5µg/ml)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.

Technical Info / Product Notes

SEPHAROSE is a trademark of GE Healthcare

companies.

UniProt ID

P18627

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-