

# LAG-3 (human) monoclonal antibody (17B4) (ATTO 647N 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- $\gamma$  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: 16

[View Online »](#)

Ordering Information

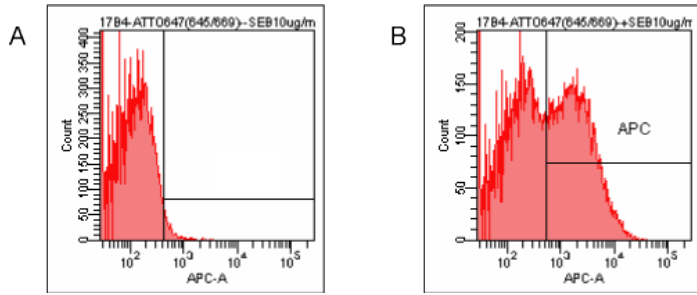
[Order Online »](#)

ALX-804-806TS0

ATTO 647N - 100 tests

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Detection of endogenous human LAG-3 by FACS analysis using LAG-3 (human), mAb (17B4) (ATTO 647) (Prod. No. ALX-804-806TS)**Method:** Human PBMC were stimulated (B) or not (A) with 1µg/ml of superantigen SEB. After 2 days PBMC were stained with 10µg/ml (1µg/0.5×10<sup>6</sup> cells) of LAG-3 (human), mAb (17B4) (ATTO 647) (Prod. No. ALX-804-806TS) and analyzed by flow cytometry.

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months after receipt when stored as recommended.
<b>Handling</b>	Avoid vigorous centrifugation and vortexing. Do not freeze. Protect from light.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Lymphocyte activation gene-3, FDC protein, CD223
<b>Application</b>	Flow Cytometry, ICC
<b>Clone</b>	17B4
<b>Formulation</b>	Liquid. In PBS containing 50mM Tris, 1% BSA and 0.02% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Synthetic peptide corresponding to 30 aa (GPPAAAPGHPLAPGPHPAAPSSWGPRPRRY) from the first N-terminal D1 domain of human LAG-3 (lymphocyte activation gene-3).
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Purified from concentrated hybridoma tissue culture supernatant by Protein A Sepharose™ CL-4B Flow Fast Column.
<b>Recommendation Dilutions/Conditions</b>	Flow Cytometry (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Source</b>	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**

**1 test means:** 1µl of MAb is used to stain 500'000 cells in a sample volume of 50µl.

ATTO-fluorescent antibodies show increased photostability, outstanding brightness and intense signals. ATTO dyes are thermally stable, resistant to environmental changes and show no significant isomerization. ATTO 647N shows red fluorescence ( $\lambda_{abs}$  (max): 645nm,  $\lambda_{em}$  (max): 669nm,  $\epsilon_{max}$ : 120'000).

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



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)