

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

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-ABS677-0100 | 100 μ g |
|-----------------|-------------|

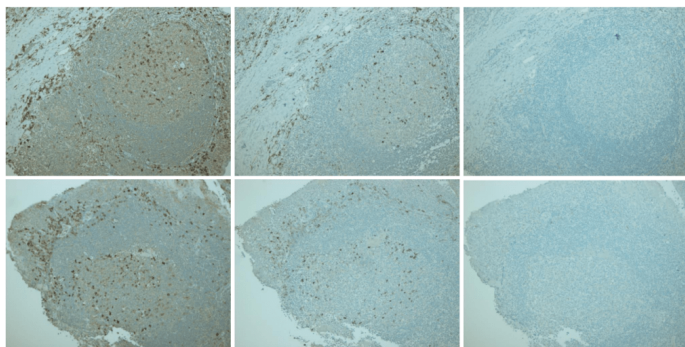
Manuals, SDS & CofA

[View Online »](#)

LAG-3 1:50

LAG-3 1:150

NO LAG-3

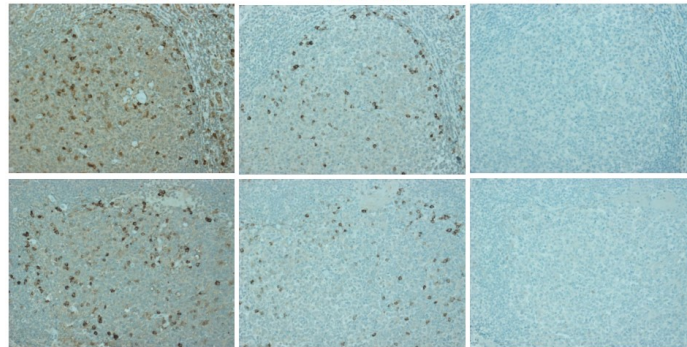


IHC of LAG-3 (human) recombinant monoclonal antibody (L4-PL33) (Prod. No. ENZ-ABS677) in human tonsil tissue (10x).

LAG-3 1:50

LAG-3 1:150

NO LAG-3

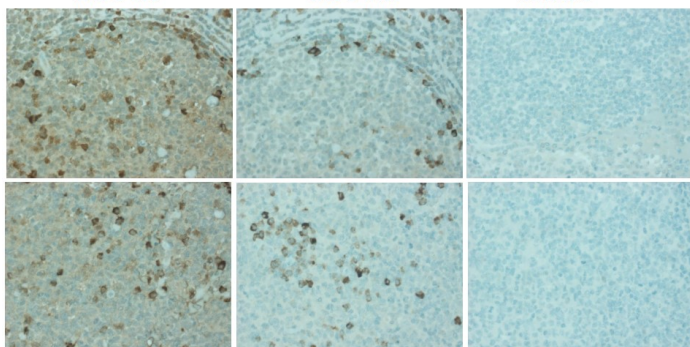


IHC of LAG-3 (human) recombinant monoclonal antibody (L4-PL33) (Prod. No. ENZ-ABS677) in human tonsil tissue (20x).

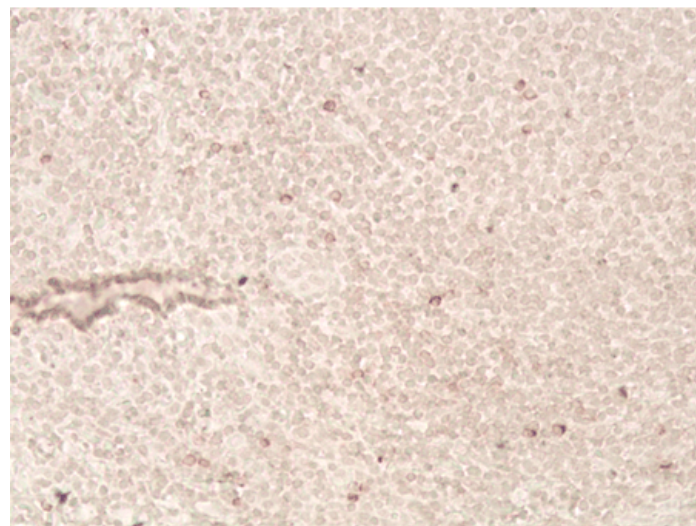
LAG-3 1:50

LAG-3 1:150

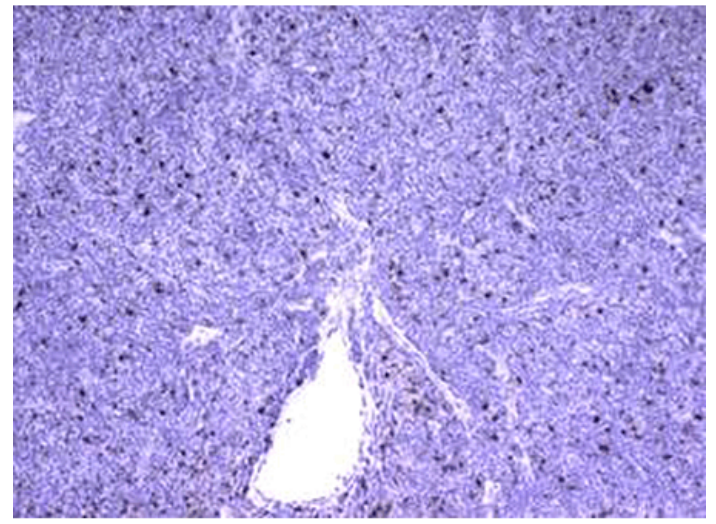
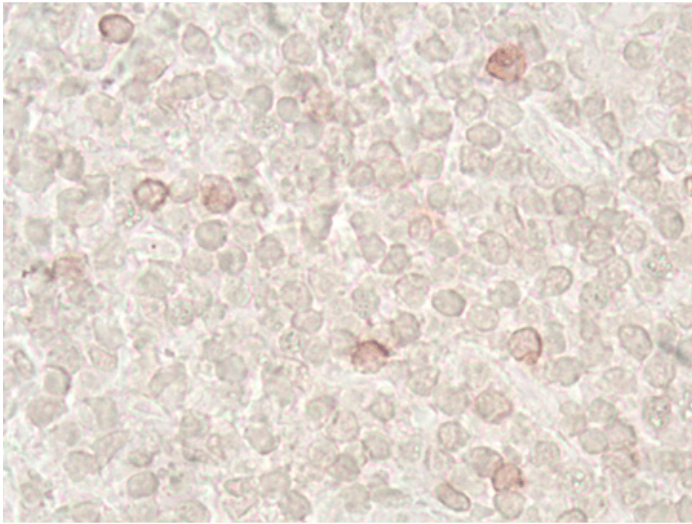
NO LAG-3



IHC of LAG-3 (human) recombinant monoclonal antibody (L4-PL33) (Prod. No. ENZ-ABS677) in human tonsil tissue (40x).

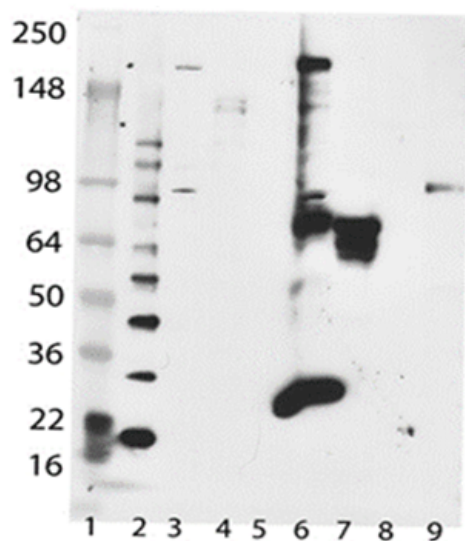


Immunohistochemistry analysis of human tonsil tissue (20x) using LAG-3 (human) recombinant mAb (L4-PL33), (Prod. No. ENZ-ABS677), at a dilution of 1:50.



Immunohistochemistry analysis of human tonsil tissue (60x) using LAG-3 (human) recombinant mAb (L4-PL33), (Prod. No. ENZ-ABS677), at a dilution of 1:50

IHC performed on human tonsil tissue using LAG-3 (human) recombinant mAb (L4-PL33) (Prod. No. ENZ-ABS677), at a dilution of 1:50.



Western blot analysis of LAG-3 recombinant mAb (L4-PL33) (Prod. No. ENZ-ABS677) using different cell lysates and purified LAG-3 protein: Lane 1 and 2: MW markers, Lane 3: Jurkat, Lane 4: HeLa, Lane 5: Empty, Lane 6: THP-2, Lane 7: HDML2, Lane 8: Empty, Lane 9: LAG-3 (human):Fc (human), (recombinant) (Prod. No. ALX-522-078).

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 containing 10% glycerol and 0.02% sodium azide. |
| 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 | IgG1κ |
| Purity Detail | Purified. |
| Recommendation Dilutions/Conditions | Immunohistochemistry (frozen sections, 1:150)Western Blot (5 µg/ml)Immunoprecipitation (10 µg/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 . |



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)