

LAG-3 (human) monoclonal antibody (11E3) (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: 4

[View Online »](#)

Ordering Information

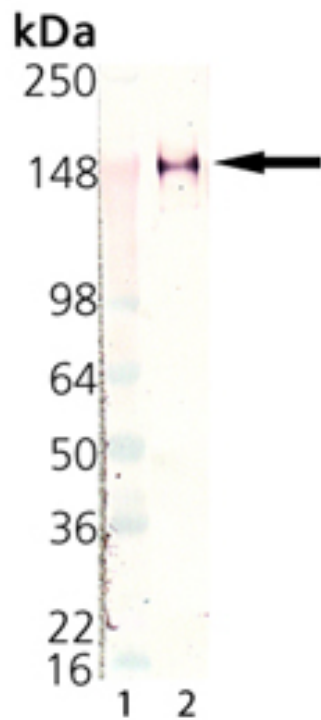
[Order Online »](#)

ALX-804-805PF-C100

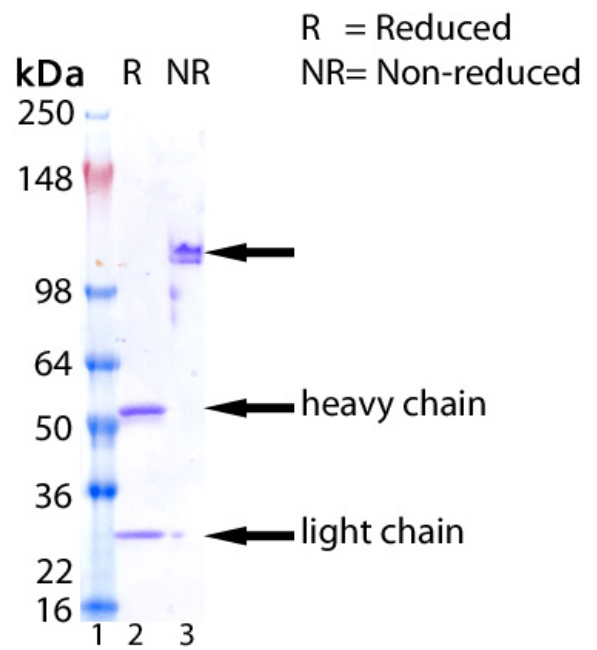
100 μ g

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of LAG-3 mAb (11E3) (Prod. No. ALX-804-805PF): Lane 1: MW marker, Lane 2: LAG-3 (human):Fc (human), (recombinant) (Prod. No. ALX-522-078).



SDS-PAGE analysis of ALX-804-805PF: Lane 1: MW Marker, Lane 2 and Lane 3: 2 μ g of LAG-3 (human) monoclonal antibody (11E3) (preservative free), resolved under reduced and non-reduced condition.

Handling & Storage

| | |
|--------------------|---------------------------|
| 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, FUNC, ICC, IHC (FS), IP, WB |
| Application Notes | Functional Application: Inhibition. |
| Clone | 11E3 |
| Formulation | Liquid. In PBS containing 10% glycerol. |
| Host | Mouse |
| Immunogen | Recombinant human LAG-3 (lymphocyte activation gene-3). |
| Isotype | IgG1 |
| Recommendation Dilutions/Conditions | ELISA (capture)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 the hybridoma culture supernatant by Protein G Fast Flow Column. |
| Species Reactivity | Human, Monkey |
| Specificity | Recognizes 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



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)