## LAG-3 (human) monoclonal antibody (11E3)

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** 

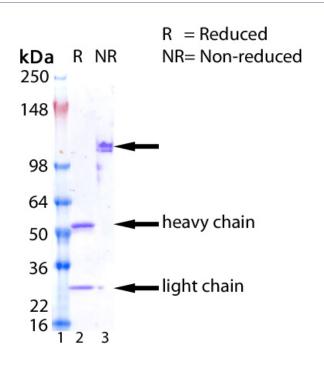
**Order Online** »

ALX-804-805-C100

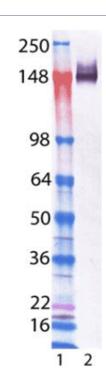
100µg

Manuals, SDS & CofA

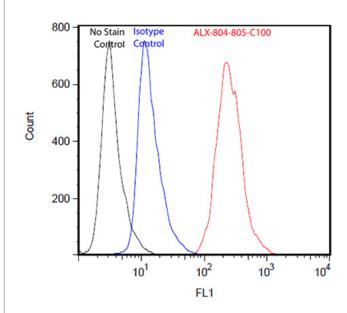
**View Online »** 



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



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



Flow cytometry analysis of 10^6 MCF-7 cells stained using LAG-3 (human) monoclonal antibody (11E3), (Prod. No. ALX-804-805), at a concentration of 50µg/mL.

## **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 NameLymphocyte activation gene-3, FDC protein, CD223

**Application** ELISA, Flow Cytometry, FUNC, ICC, IHC (FS), IP, WB

Clone 11E3

Formulation Liquid. In PBS containing 10% glycerol and 0.02% sodium

azide.

**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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com