

TL1A (human) monoclonal antibody (L4G9)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

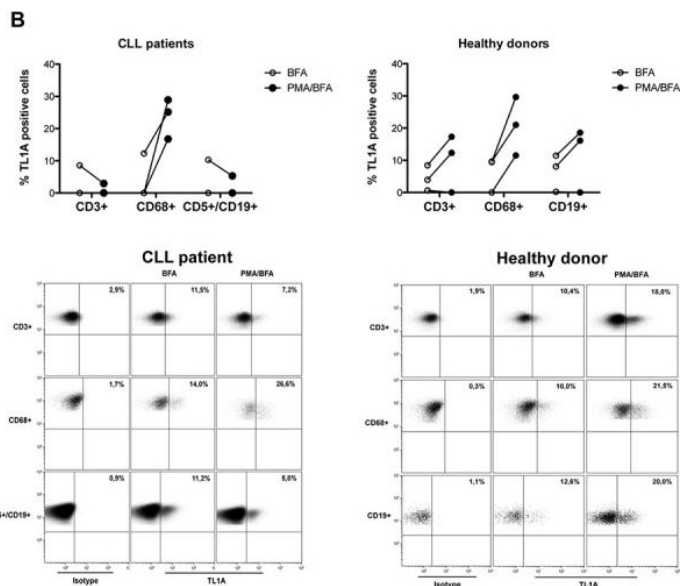
Ordering Information

[Order Online »](#)

ALX-804-832-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



TL1A expression in PBMC cell subsets from CLL patients and healthy donors. A. Quantitative real-time PCR analysis of TL1A mRNA expression in FACS-isolated cell subsets (CD3+, CD14+, and CD19+) from CLL (n = 6) and healthy donors (n = 3). Gene expression was normalized vs β -actin expression and represented as relative expression of TL1A mRNA levels with respect to mean TL1A mRNA across samples (set as 1). Lines represent median values. B. TL1A protein expression analysis in gated-cell subsets (CD3+, CD68+, CD5+/CD19+, CD19+) of PBMCs from CLL patients (n = 3) and healthy donors (n = 3). PBMCs were unstimulated or stimulated with PMA/ionomycin (PMA) in presence of brefeldin A (BFA). Dot plots are representative examples of TL1A flow cytometry analysis in gated-cell subsets (CD3+, CD68+, CD5+/CD19+, CD19+) of PBMCs from CLL patient and healthy donor. Number in bold indicate the percentage of TL1A-positive cells.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Expression and function of the TL1A/DR3 axis in chronic lymphocytic leukemia. *Oncotarget* (2015)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TNFSF 15, VEGI
Application	ELISA, Flow Cytometry
Clone	L4G9
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human TL1A extracellular domain.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human

UniProt ID

O95150



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Worry-free Guarantee

This antibody is covered by our **Worry-Free Guarantee**.