

TL1A (soluble) (mouse), (recombinant)

TL1A (TNFSF15), a ligand belonging to the tumor necrosis factor (TNF) family, is expressed predominantly by endothelial cells and monocytes. TL1a is upregulated by the proinflammatory cytokines TNF and IL-1 and also by immune complexes (IC). TL1 functions in T cell costimulation and Th1 polarization. On activated T cells, TL1A functions specifically via its surface-bound receptor DR3, (a member of the death-domain containing TNF receptor family) to promote cell survival and secretion of proinflammatory cytokines. The secreted decoy receptor 3 (DcR3), a soluble protein of the tumor necrosis factor receptor (TNFR) superfamily blocks the action of TL1A. TL1A can be released to circulate as a homotrimeric soluble form. Activation of DR3 by TL1A induced the formation of a signaling complex containing TRADD, TRAF2, and RIP and activated the NF- κ B and the ERK, JNK, and p38 mitogen-activated protein kinase pathways. The TL1A/DR3 pathway plays an important role in Th1-mediated intestinal diseases, such as Crohn's disease.

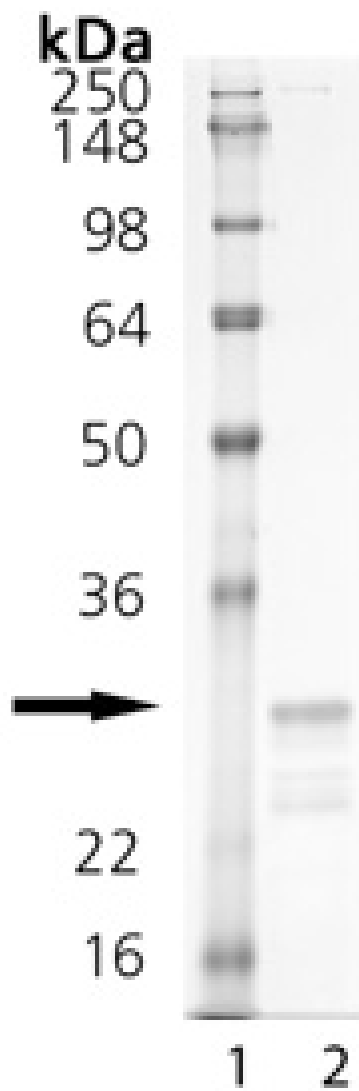
Ordering Information

[Order Online »](#)

ALX-522-105-C010	10 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of TL1A (soluble) (mouse), (recombinant), Lane 1: MW Marker, Lane 2: 1 μ g mouse TL1A. Banding observed beneath primary band is immunoreactive with anti-FLAG antibody.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TNFSF 15, VEGI
Application Notes	ELISA: binds to DcR3. Flow Cytometry: use together with Enhancer for Ligands (Prod. No. ALX-804-034) to detect human and mouse DR3 and DcR3.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~25kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of mouse TL1A (aa 94-252) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	Q8K3Y8



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)