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-kB 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

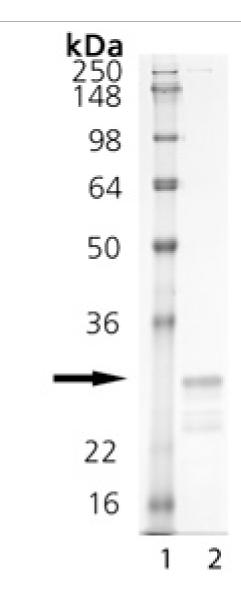
Order Online »

ALX-522-105-C010

10µg

Manuals, SDS & CofA

View Online »



SDS-PAGE analysis of TL1A (soluble) (mouse), (recombinant), Lane 1: MW Marker, Lane 2: 1µg 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)

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com