

TNC:TWEAK (soluble) (mouse), (recombinant)

Some TNF ligands (such as TWEAK (mouse) or FasL (mouse)) produced as soluble forms do not assemble into stable trimers and show low receptor binding. Addition of the tenascin-C oligomerization domain (TNC) (from chicken) at the N-terminus of recombinant mouse soluble TWEAK stabilizes the TNF homology domain (THD) of rmsTWEAK into a trimer and increases binding to its respective TNF receptor. Note: For some TNF ligands, such as FasL (mouse), activated by further oligomerization (> hexamer), the presence of TNC is not sufficient and ligand crosslinking using the "Enhancer for Ligands" (Prod. No. ALX-804-034) is required.

Ordering Information

[Order Online »](#)

ALX-522-138-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

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.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Apo-3L, TNFSF 12
Application Notes	ELISA: binds to mouse Fn14.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Associates of CAPE COD Inc.).
Formulation	Lyophilized. Contains PBS.
MW	~30kDa
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 TWEAK (aa 94-249) is fused at the N-terminus to tenascin-C (TNC) (chicken) (aa 110-139) and a FLAG [®] -tag.
Specificity	Binds to Fn14 (mouse):Fc (human) (rec.) (Prod. No. ALX-522-036)

Technical Info / Product FLAG is a registered trademark of Sigma-Aldrich Co.

Notes



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

054907

European Sales Office
ENZO LIFE SCIENCES
(EU) 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