

TNF-R1 (human):Fc (human), (recombinant)

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-013-C050	50µg
------------------	------

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.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	p55, CD120a, TNF-R 60kDa, TNFRSF 1A
Application Notes	ELISA: binds to TNF- α .
Concentration	1mg/ml after reconstitution.
Formulation	Lyophilized. Contains PBS.
MW	~60kDa (reducing SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50 μ l sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of human TNF-R1 (aa 1-211) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse TNF- α (tumor necrosis factor- α).
Technical Info / Product Notes	Previous data has shown that TNF-R1 receptor binds to TNF-R1 and triggers various biological effects (including cell death). rhTNF-R1:Fc fusion protein blocks TNF- α -mediated lysis with a inhibitory activity in a concentration range of 10-50ng/ml using Wehi cells.
UniProt ID	P19438

Last modified: May 29, 2024



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)