

Fc (human):TNF- α (soluble) (human), (recombinant)

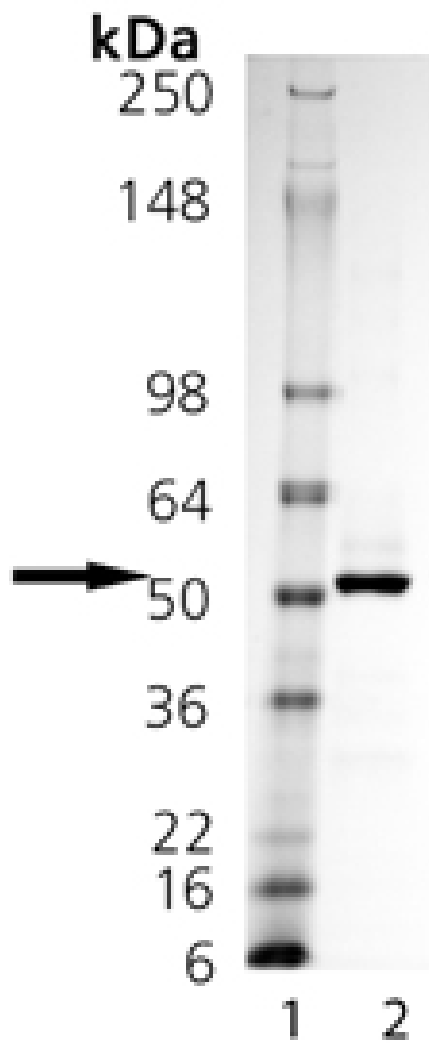
Ordering Information

[Order Online »](#)

ALX-522-087-C010	10 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of Fc (human):TNF- α (soluble) (human), (recombinant): Lane 1: MW Marker, Lane 2: 1 μ g TNF- α .

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	Tumor necrosis factor α , TNFSF 2
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/ μ g purified protein (LAL test; Associates of CAPE COD Inc.).
Formulation	Lyophilized. Contains PBS.
MW	~48kDa (calculated), ~52kDa (observed)
Purity	\geq 90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute with 100 μ l sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in CHO cells. The extracellular domain of human TNF- α (aa 85-233) is fused at the N-terminus to the Fc portion of human IgG1 and a linker peptide (20 aa).
Technical Info / Product Notes	Historical data has shown that TNF α induces killing of WEHI-164 cells in a concentration range of 0.5-2ng/ml (does not require an enhancer).
UniProt ID	P01375



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)