

TNF- α (soluble) (mouse), (recombinant)

Citations: 22

[View Online »](#)

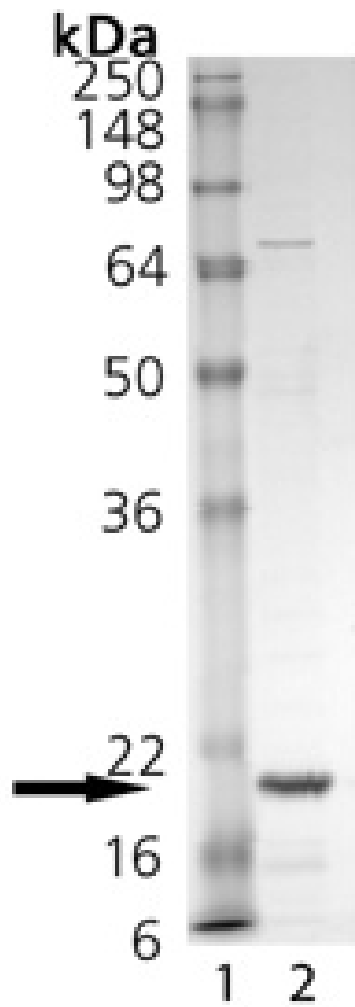
Ordering Information

[Order Online »](#)

ALX-522-009-C050	50 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis of TNF- α (soluble) (mouse), (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
Application Notes	ELISA: binds to human TNF-R1 receptor.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/ μ g purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
Gene/Protein Identifier	A002291 (UCSD Signaling Gateway ID)
MW	~20kDa (SDS-PAGE).
Purity	\geq 90% (SDS-PAGE)
Reconstitution	Reconstitute with 50 μ l sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in <i>E. coli</i> . The extracellular domain of mouse TNF- α (tumor necrosis factor- α) (aa 77-235) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Specificity	Binds to human, mouse and rat TNF-R1 and less efficiently to TNF-R2. In the presence of cross-linking enhancer (see Set Prod. No. ALX-850-061), TNF- α shows a significantly higher affinity for TNF-R2 than for TNF-R1, mimicking the characteristics of membrane-bound TNF- α .

Technical Info / Product Notes	<p>Historical lots have shown that it exerts its biological activity in a concentration range of 0.5-1ng/ml (WEHI 164 cells).</p> <p>FLAG is a registered trademark of Sigma-Aldrich Co.</p>
UniProt ID	P06804



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) 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