

# TNF- $\alpha$ (soluble) (mouse), (recombinant)

Citations: 22

[View Online »](#)

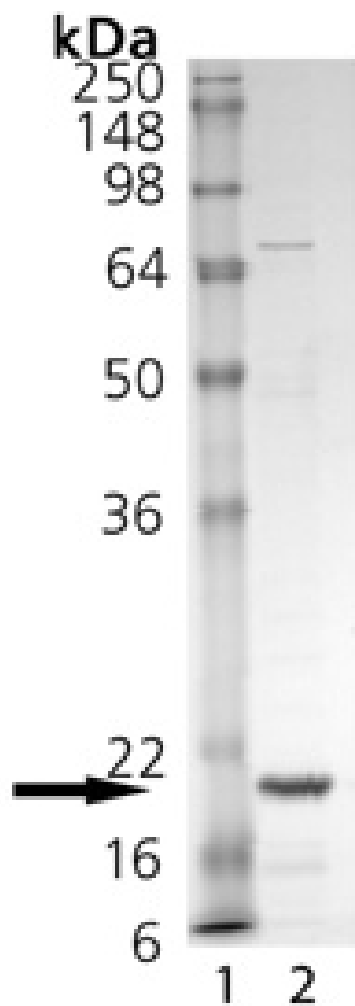
## Ordering Information

[Order Online »](#)

ALX-522-009-C050	50 $\mu$ g
------------------	------------

## Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis of TNF- $\alpha$  (soluble) (mouse), (recombinant). Lane 1: MW Marker, Lane 2: 1  $\mu$ g TNF- $\alpha$

# 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/μg 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	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 50μl 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 <sup>®</sup> -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

Historical lots have shown that it exerts its biological activity in a concentration range of 0.5-1ng/ml (WEHI 164 cells).

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

P06804

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)