

# TNF- $\alpha$ (human), (recombinant) (cell culture grade)

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-520-0020	10 $\mu$ g
--------------	------------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Lyophilized protein is stable at room temperature for a few weeks.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Tumor necrosis factor- $\alpha$ , TNFSF 2
<b>Biological Activity</b>	Induces cytolysis of mouse L929 cells in the presence of 1 $\mu$ g/ml Actinomycin D (Prod. No. BML-GR300) (ED <sub>50</sub> $\leq$ 0.05ng/ml).
<b>CAS</b>	94948-59-1
<b>Formulation</b>	Lyophilized from 0.2 $\mu$ m-filtered solution containing mannitol and trehalose.
<b>MI</b>	<b>14:</b> 9812
<b>MW</b>	~17kDa.
<b>Purity</b>	$\geq$ 97% (SDS-PAGE)
<b>Reconstitution</b>	A quick spin is recommended followed by reconstitution in distilled, sterile water to a concentration of 0.1-1mg/ml. For further dilution, add a carrier protein to avoid loss of biological activity.
<b>Source</b>	Produced in yeast.
<b>Specific Activity</b>	$\geq 2 \times 10^7$ U/mg protein. Determined by standard cell cytotoxicity assay (N. Matthes & M.L. Neale; Lymphokines and Interferons, A Practical Approach; M.J. Clemens et al., eds.; IRL Press, p. 221).
<b>UniProt ID</b>	P01375

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)