

TNF- α (human), (recombinant) (cell culture grade)

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-520-002-C010	10 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Tumor necrosis factor- α , TNFSF 2
Biological Activity	Induces cytolysis of mouse L929 cells in the presence of 1 μ g/ml Actinomycin D (Prod. No. BML-GR300) (ED ₅₀ \leq 0.05ng/ml).
CAS	94948-59-1
Formulation	Lyophilized from 0.2 μ m-filtered solution containing mannitol and trehalose.
MI	14: 9812
MW	~17kDa.
Purity	\geq 97% (SDS-PAGE)
Reconstitution	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.
Source	Produced in yeast.
Specific Activity	$\geq 2 \times 10^7$ 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).
UniProt ID	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)