

TNF-R2 (human):Fc (human), (recombinant)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-0140	50µg
--------------	------

Manuals, SDS & CofA

[View Online »](#)

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	p75, CD120b, TNF-R 80kDa, TNFRSF 1B
Application Notes	ELISA: binds to TNF- α . Detection of membrane-bound TNF- α in combination with PAb to human IgG1.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/ μ g purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~75kDa (reducing SDS-PAGE).
Purity	\geq 95% (SDS-PAGE)
Reconstitution	Reconstitute with 50 μ l sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of human TNF-R2 (aa 1-257) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse TNF- α (tumor necrosis factor- α).
UniProt ID	P20333

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)