

# TNF-R1 (human):Fc (human), (recombinant)

Citations: 8

[View Online »](#)

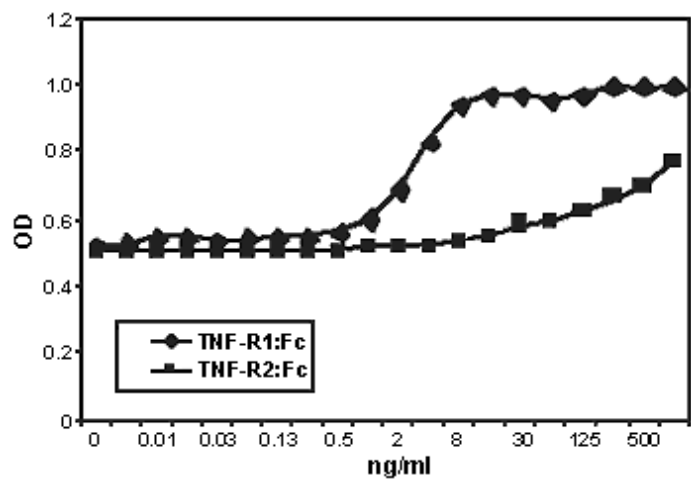
## Ordering Information

[Order Online »](#)

ALX-522-013-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Inhibition of TNF-mediated lysis of WEHI 164 target cells (rhTNF-a, Prod. No. [ALX-522-008](#)). rhTNF-R1:Fc exerts its inhibitory activity in a concentration range of 10-50ng/ml.””,TNF-R1 (human):Fc (human)

## 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	p55, CD120a, TNF-R 60kDa, TNFRSF 1A
Application Notes	ELISA: binds to TNF- $\alpha$ .
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/ $\mu$ g purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~60kDa (reducing SDS-PAGE).
Purity	$\geq$ 95% (SDS-PAGE)
Reconstitution	Reconstitute with 50 $\mu$ l sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of human TNF-R1 (aa 1-211) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse TNF- $\alpha$ (tumor necrosis factor- $\alpha$ ).
Technical Info / Product Notes	Previous data has shown that TNF-R1 receptor binds to TNF-R1 and triggers various biological effects (including cell death). rhTNF-R1:Fc fusion protein blocks TNF- $\alpha$ -mediated lysis with a inhibitory activity in a concentration range of 10-50ng/ml using Wehi cells.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

P19458

European Sales Office  
ENZO LIFE SCIENCES  
(UK) 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-fr@enzolifesciences.com](mailto:info-fr@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)