

# TNF receptor 1

## polyclonal antibody

Tumor necrosis factor alpha (TNF- $\alpha$ ), also known as cachectin, is a 17.5 kDa 157 amino acid member of the TNF superfamily of cytokines that is a potent lymphoid factor with effects on a wide range of target cells. Active TNF $\alpha$  is produced in both soluble and membrane-anchored trimers by macrophages, NK cells, and T- and B-lymphocytes. TNF exerts proinflammatory signals via binding and inducing trimerization of TNF-receptor1 (TNF-R1) expressed on most normal and transformed cells, or to TNF-receptor 2 (TNF-R2), expressed on endothelial and most immune cells. TNF signaling regulates hematopoiesis, differentiation, endothelial cell activation, apoptosis, lipid metabolism, tumor progression, and immune surveillance, and dysregulation of TNF or its receptors is implicated in numerous disease states including cancer, osteoporosis, autoimmune disease, diabetes, and atherosclerosis.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 8

[View Online »](#)

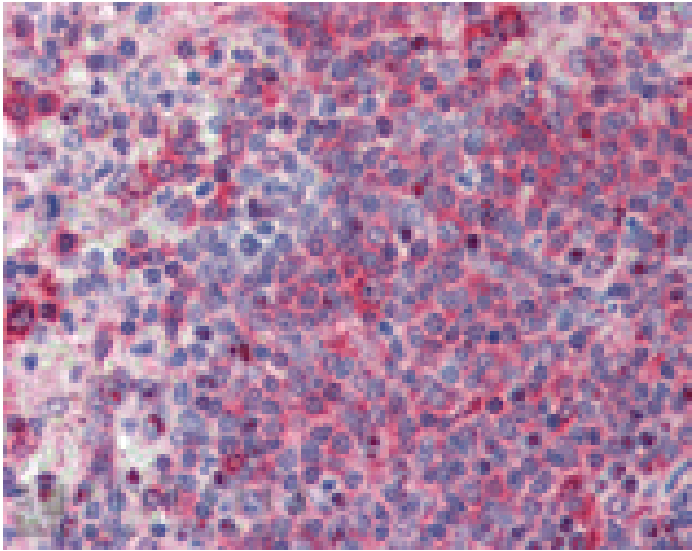
### Ordering Information

[Order Online »](#)

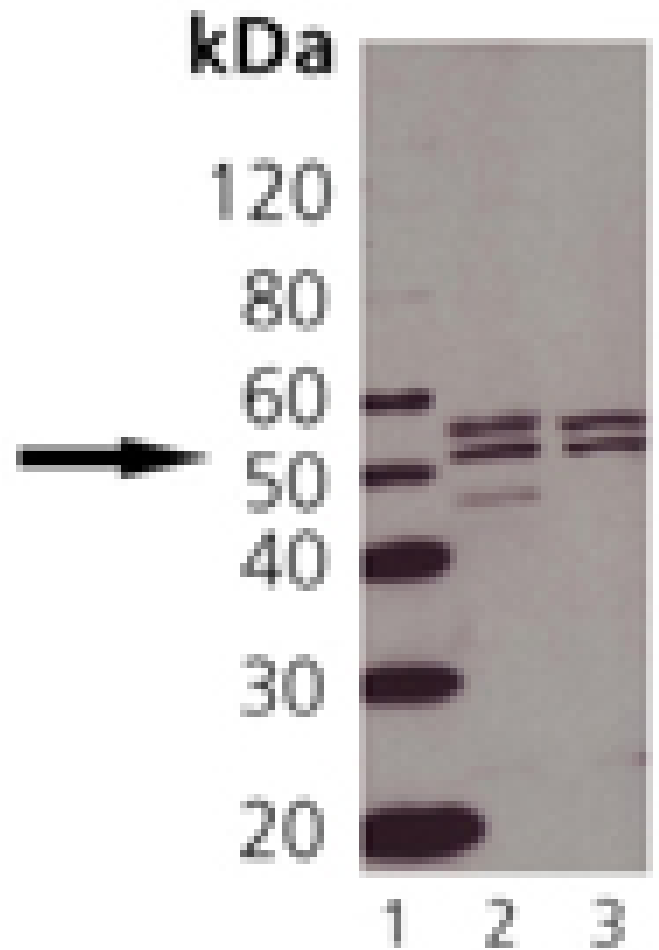
ADI-CSA-815-D	50 $\mu$ g
ADI-CSA-815-F	200 $\mu$ g

### Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human spleen tissue stained with TNF receptor 1 pAb at 10µg/ml.



Western blot analysis of TNF-R1: Lane 1: MW marker, Lane 2: Hela (HS), Lane 3: Jurkat

# Handling & Storage

Long Term Storage                    -20°C

Shipping                                Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Tumor necrosis factor receptor 1, TNF-R1
Application	IHC (PS), IP, WB
Application Notes	Detects a band of 55kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	X59238
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of mouse TNF-R1. The sequence is completely conserved in rat and human.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (12.5µg/ml)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Bovine, Dog, Drosophila, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus, Yeast
UniProt ID	P25118
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



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)