

# TNF- $\alpha$ Recombinant monoclonal antibody (cA2) (Human IgG1 $\kappa$ )

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

## Ordering Information

[Order Online »](#)

ENZ-ABS562-0200	200 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	TNFSF 2, Tumor necrosis factor- $\alpha$ , Cachectin, Tumor necrosis factor ligand superfamily member 2
Application	ELISA, Flow Cytometry, in vivo Assay
Clone	cA2 (Infliximab)
Crossreactivity	Does not cross-react with TNF- $\beta$ .
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Human
Immunogen	Purified recombinant human TNF (rTNF).
Isotype	IgG1 $\kappa$
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	Binds to soluble and transmembrane forms of TNF- $\alpha$ and neutralizes its biological activity.
Technical Info / Product Notes	This is a chimeric antibody created for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P01375

Worry-free Guarantee

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



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)