

# CD4 Recombinant monoclonal antibody (CE9.1) (Human IgG4 $\lambda$ )

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

## Ordering Information

[Order Online »](#)

ENZ-ABS440-0200	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	T-cell differentiation antigen, Ly-4, L3T4, T-cell surface antigen T4/Leu-3
Application	ELISA, Flow Cytometry, FUNC, IHC
Clone	CE9.1 (Clenoliximab)
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Human
Immunogen	Domain 1 of human CD4.
Isotype	IgG4λ
Purity Detail	Protein A affinity purified.
Species Reactivity	Chimpanzee, Human
Specificity	The antibody binds to human CD4 with a Kd of 1.0 nM and an ED50 of 0.2 µg/ml.
Technical Info / Product Notes	This reformatted human antibody was made using the variable domain sequences of the original Human IgG4 (primatized) format, for improved compatibility with existing reagents, assays and techniques.



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)

**Worry-free Guarantee**

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

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.