

CD4 Recombinant monoclonal antibody (CE9.1) (Rabbit IgG λ)

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

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	Rabbit
Immunogen	Domain 1 of human CD4.
Isotype	IgG λ
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 μ g/ml.
Technical Info / Product Notes	This chimeric rabbit antibody was made using the variable domain sequences of the original Human IgG4 (primatized) format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P01730

Worry-free Guarantee

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

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)