## CD4 Recombinant monoclonal antibody (CE9.1) (Rabbit IgGλ)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS439-0200

200µg

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\(\lambda\)

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  $\mu$ 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.

P01730

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



eu@enzolifesciences.com