CD4 Recombinant monoclonal antibody (MT310) (Rabbit IgGк)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS443-0200

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 Flow Cytometry, IF, IHC (PS), IP, WB

Clone MT310

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Isotype IgGĸ

Purity Detail Protein A affinity purified.

Species Reactivity Human

Specificity Recognizes the extracellular domain D1 of CD4 – the

same site as the HIV-1 gp120-binding site.

Technical Info / Product NotesThis chimeric rabbit antibody was made using the variable

domain sequences of the original murine IgG1 format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID P01730

Last modified: May 29, 2024



info-

eu@enzolifesciences.com