CD4 Recombinant monoclonal antibody (YNB46.1.8) (Rabbit IgGk)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS430-0200

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, IHC (FS)

Clone YNB46.1.8 (Campath-9H)

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen Rat T-cell line (NB2) transfected with the human CD4.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

Specificity Recognizes the 55kDa human CD4 cell surface antigen,

expressed by helper T lymphocytes and by monocytes.

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

domain sequences of the original rat IgG1 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