CD3ε Recombinant monoclonal antibody (145-2C11) (Rabbit IgGκ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS424-0200

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 surface glycoprotein CD3 ε chain, T-cell surface

antigen T3/Leu-4 ε chain, CD3e

Application Flow Cytometry, FUNC, IP, WB

Application Notes Functional Application: Activation

Clone 145-2C11

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen H-2Kb-specific murine cytotoxic T-lymphocyte (CTL) clone.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Mouse, Rat

Specificity Recognizes a 25-kDa protein component (CD3e, originally

called T3-e) of the antigen specific T-cell receptor.

Technical Info / Product Notes

This chimeric rabbit antibody was made using the variable domain sequences of the original Hamster IgG1 format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P22646

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-