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

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS425-0200

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.

lgGk

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.

UniProt ID P22646

Last modified: May 29, 2024



info-

eu@enzolifesciences.com