CD3ε Recombinant monoclonal antibody (12F6) (Rabbit IgGκ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS622-0200

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 epsilon chain, CD3e

Application Flow Cytometry

Clone 12F6

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen Human CD3.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

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

domain sequences of the original Mouse IgG2a format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID P07766

Last modified: May 29, 2024



info-

eu@enzolifesciences.com