## CD3ε Recombinant monoclonal antibody (UCHT1) (Mouse IgG1κ)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS429-0200

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, IHC

Clone UCHT1

**Formulation** Liquid. In PBS containing 0.02% Proclin 300.

**Host** Mouse

**Immunogen** Human infant thymocytes.

IgG1ĸ

Purity Detail Protein A affinity purified.

Species Reactivity Human

**Specificity** Recognizes epitopes on CD3-ε found in heterodimers with

CD3- $\delta$  and with CD3- $\gamma$  of  $\alpha/\beta$  T cell receptor (TCR).

UniProt ID P07766

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com