## CD3ε Recombinant monoclonal antibody (YTH 12.5) (Rabbit IgGλ)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS624-0200

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

Clone YTH 12.5

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

**Host** Rabbit

Immunogen Human CD3.

lgGλ

Purity Detail Protein A affinity purified.

Species Reactivity Human

**Technical Info / Product Notes**This chimeric rabbit antibody was made using the variable

domain sequences of the original rat IgG2b format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID P07766

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com