## CD8 β Recombinant monoclonal antibody (YTS 156.7.7) (Rat IgG2bκ)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS631-0200

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 CD8 beta chain, Lymphocyte

antigen 3, Ly-3

**Application** Flow Cytometry, IHC

**Clone** YTS 156.7.7

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

**Host** Rat

Immunogen Mouse CD8 (lyt-3).

Isotype IgG2bk

Purity Detail Protein A affinity purified.

Species Reactivity Mouse

**Specificity** The antibody recognizes the B epitope of the mouse Lyt-3

(CD8) antigen. Has been used in CD8 subset depletion in

in vivo and flow cytometry studies.

UniProt ID P10300

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com