CD8 β Recombinant monoclonal antibody (YTS 156.7.7) (Rabbit IgGκ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS632-0200

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-4

Application Flow Cytometry, IHC

Clone YTS 156.7.7

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen Mouse CD8 (lyt-3).

Isotype IgGk

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.

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

domain sequences of the original format, for improved

compatibility with existing reagents, assays and

techniques.

P10300

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com