

# CD8 $\beta$ Recombinant monoclonal antibody (YTS 156.7.7) (Rat IgG2b $\kappa$ )

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

ENZ-ABS631-0200	200 $\mu$ g
-----------------	-------------

## 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 (Iyt-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 <i>in vivo</i> and flow cytometry studies.
UniProt ID	P10300

### Worry-free Guarantee



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

### This antibody is covered by our **Worry-Free Guarantee.**