

# CD8 $\alpha$ Recombinant monoclonal antibody (YTS 105.18) (Rabbit IgG $\kappa$ )

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

## Ordering Information

[Order Online »](#)

ENZ-ABS456-0200	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	CD8a, OKT8 T-cell antigen, Leu2 T-lymphocyte antigen, T8 T-cell antigen, p32, T-cell surface glycoprotein CD8 α chain, T-cell surface glycoprotein Lyt-2
Application	Flow Cytometry, FUNC, IF, IHC (FS)
Application Notes	This antibody has been reported to block MHC I dependent T cell responses <i>in vitro</i> and <i>in vivo</i> .
Clone	YTS 105.18
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	Ly-2-transfected mouse L cells.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Species Reactivity	Mouse
Specificity	Recognizes a non polymorphic epitope on the mouse CD8 α chain.

## Technical Info / Product Notes

This chimeric rabbit antibody was made using the variable domain sequences of the original rat IgG2a format, for improved compatibility with existing reagents, assays and techniques.

### UniProt ID

P01731

### Worry-free Guarantee

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

Last modified: May 29, 2024



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)