

CD8 α Recombinant monoclonal antibody (YTS 105.18) (Rabbit IgG κ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS456-0200	200 μ 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	IgGk
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](#).



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)