

CD8 α Recombinant monoclonal antibody (YTS 105.18) (Rat IgG2 κ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS457-0200	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	Rat
Immunogen	Ly-2-transfected mouse L cells.
Isotype	IgG2ak
Purity Detail	Protein A affinity purified.
Species Reactivity	Mouse
Specificity	Recognizes a non polymorphic epitope on the mouse CD8 α chain.
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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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