

CD3ε Recombinant monoclonal antibody (YTH 12.5) (Rat IgG2bλ)

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

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS623-0200 | 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 CD3 ε chain, T-cell surface antigen T3/Leu-4 epsilon chain, CD3e |
| Clone | YTH 12.5 |
| Formulation | Liquid. In PBS containing 0.02% Proclin 300. |
| Host | Rat |
| Immunogen | Human CD3. |
| Isotype | IgG2bλ |
| Purity Detail | Protein A affinity purified. |
| Species Reactivity | Human |
| UniProt ID | P07766 |
| 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-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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