

# CD3 (human) monoclonal antibody (TR66)

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

Citations: 12

[View Online »](#)

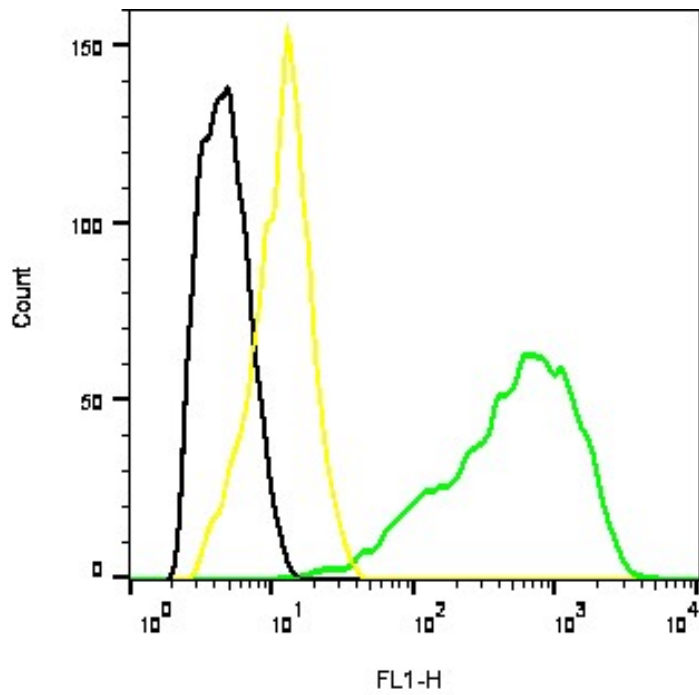
## Ordering Information

[Order Online »](#)

ALX-804-822-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of  $10^6$  Jurkat cells using the isotype control of IgG1 (Prod. No. ADI-SAB-600) compared with CD3 (human) mAb (Prod. No. ALX-804-822). Cells were preincubated with unconjugated antibody and then stained with conjugated secondary antibody (Invitrogen, Cat#: A11001).

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	For long term storage, aliquot and freeze at -20°C. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	Flow Cytometry, FUNC, ICC, IP, WB
Application Notes	Detects endogenous protein by FC, ICC and IP.
Clone	TR66
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Cloned human helper T cell line.
Isotype	IgG1
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human
Specificity	Recognizes the $\epsilon$ -chain of CD3.
UniProt ID	P09693 ( $\gamma$ chain), P04234 ( $\delta$ chain), P07766 ( $\epsilon$ chain), P20963 ( $\zeta$ chain)

### 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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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

Last modified: May 29, 2024