

# CD3 $\epsilon$ Recombinant monoclonal antibody (145-2C11) (Mouse IgG1 $\kappa$ )

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

## Ordering Information

[Order Online »](#)

ENZ-ABS426-0200	200 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	T-cell surface glycoprotein CD3 ε chain, T-cell surface antigen T3/Leu-4 ε chain, CD3e
<b>Application</b>	Flow Cytometry, FUNC, IP, WB
<b>Application Notes</b>	Functional Application: Activation
<b>Clone</b>	145-2C11
<b>Formulation</b>	Liquid. In PBS containing 0.02% Proclin 300.
<b>Host</b>	Rabbit
<b>Immunogen</b>	H-2Kb-specific murine cytotoxic T-lymphocyte (CTL) clone.
<b>Isotype</b>	IgG1κ
<b>Purity Detail</b>	Protein A affinity purified.
<b>Species Reactivity</b>	Mouse, Rat
<b>Specificity</b>	Recognizes a 25-kDa protein component (CD3e, originally called T3-e) of the antigen specific T-cell receptor.

## Technical Info / Product Notes

This chimeric mouse antibody was made using the variable domain sequences of the original Hamster IgG1 format, for improved compatibility with existing reagents, assays and techniques.

## UniProt ID

P22646

## 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)