

CD3ε Recombinant monoclonal antibody (145-2C11) (Rabbit IgGκ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS424-0200	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 ε chain, CD3e
Application	Flow Cytometry, FUNC, IP, WB
Application Notes	Functional Application: Activation
Clone	145-2C11
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	H-2Kb-specific murine cytotoxic T-lymphocyte (CTL) clone.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Species Reactivity	Mouse, Rat
Specificity	Recognizes a 25-kDa protein component (CD3e, originally called T3-e) of the antigen specific T-cell receptor.
Technical Info / Product Notes	This chimeric rabbit 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](#).



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)