

CD3ε Recombinant monoclonal antibody (12F6) (Mouse IgG2ak)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS621-0200	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
Application	Flow Cytometry
Clone	12F6
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Human CD3.
Isotype	IgG2ak
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-be@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-fr@enzolifesciences.com

France
Phone: +33 472 440 655

info-de@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