

# CD3ε recombinant monoclonal antibody (RM344)

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

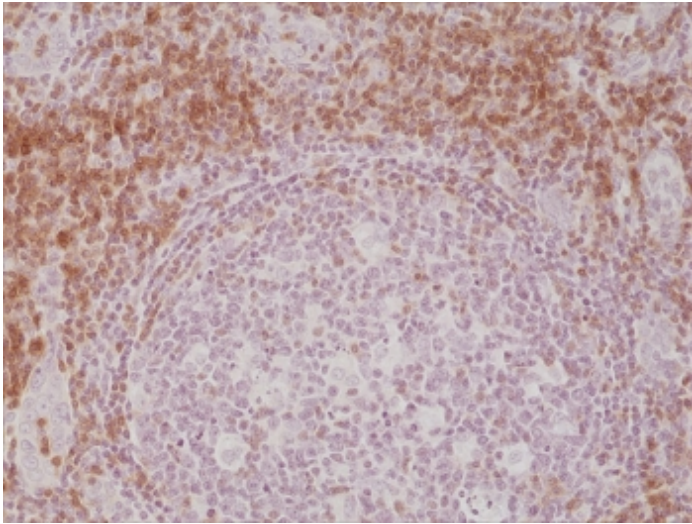
## Ordering Information

[Order Online »](#)

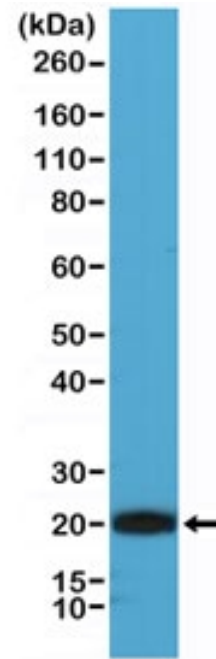
ENZ-ABS878-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



IHC staining of FFPE human tonsil tissue with recombinant CD3 epsilon antibody at 1:200.



Western blot testing of human Jurkat cell lysate with recombinant CD3 epsilon antibody at 1:1000. Predicted molecular weight ~23 kDa.

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      T-cell surface antigen T3/Leu-4 epsilon chain

Application      IHC (PS), WB

Clone      RM344

Formulation      Liquid. In PBS containing 50% glycerol, 1% BSA and 0.09% sodium azide.

Host      Rabbit

Immunogen      Peptide corresponding to the C-terminus of human CD3ε chain.

Isotype      IgG

Purity Detail      Protein A purified from animal origin-free supernatant.

**Recommendation**  
**Dilutions/Conditions**      Immunohistochemistry (FFPE) (1:100-1:200)Western blot (1:1,000-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity      Human

UniProt ID      P07766

**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-  
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)