## CD3ε recombinant monoclonal antibody (RM344)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

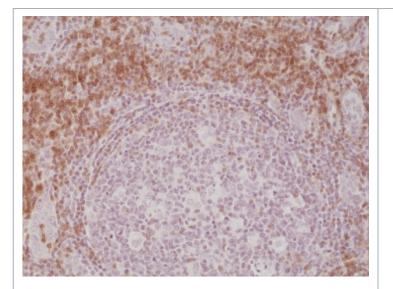
Order Online »

**ENZ-ABS878-0100** 

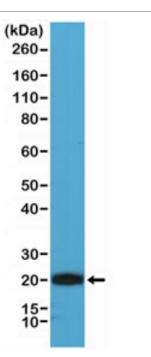
100µl

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

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com