

# CD5 Recombinant monoclonal antibody (RM314)

T-cell surface glycoprotein CD5 may act as a receptor in regulating T-cell proliferation.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS672-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** T-cell surface glycoprotein CD5, Lymphocyte antigen T1/Leu-1

**Application** IHC (PS), WB

**Clone** RM314

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

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the C-terminus of human CD5.

**Isotype** IgG

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (FFPE): 1:500-1:1,000 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.

**Source** Purified from animal origin-free supernatant.

**Species Reactivity** Human

**UniProt ID** P06127

**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)