

# CD40 (human) monoclonal antibody (mAb 89) (biotin conjugate)

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

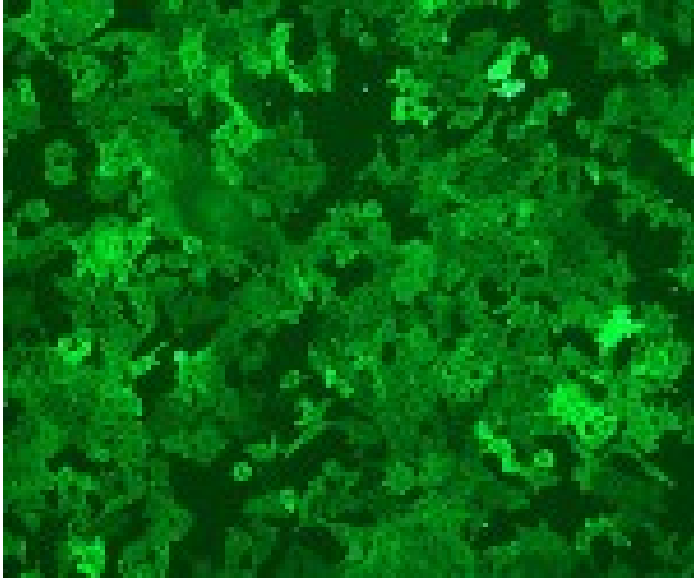
## Ordering Information

[Order Online »](#)

ENZ-ABS146-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of CD40-transfected L cells stained with CD40 (human) mAb, clone mAb89.

## Handling & Storage

**Short Term Storage** -20°C

**Long Term Storage** -80°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Tumor necrosis factor receptor superfamily member 5, TNFRSF 5, B-cell surface antigen CD40, Bp50, CD40L receptor, CDw40

**Application** Flow Cytometry, IHC (PS)

**Clone** mAb 89

**Formulation** Liquid. In PBS containing 50% glycerol.

**Host** Mouse

**Immunogen** Anti-μ activated tonsillar human B cells.

**Isotype** IgG1

**Purity Detail** Purified using QMA Hyper D ion exchange chromatography

**Species Reactivity** Human

**UniProt ID** P25942

**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)