

# MAX recombinant monoclonal antibody (103)

MAX (Myc-associated factor X) is a transcription factor that dimerizes with MYC and other proteins to regulate gene expression involved in cell growth, apoptosis, and differentiation. It plays a crucial role in oncogenesis and is studied in cancer research for its interactions with the MYC signaling pathway.

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

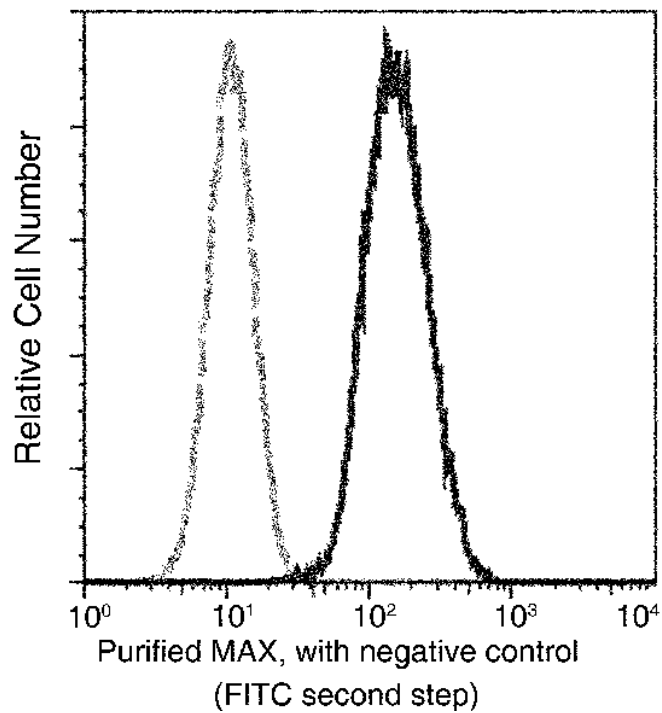
## Ordering Information

[Order Online »](#)

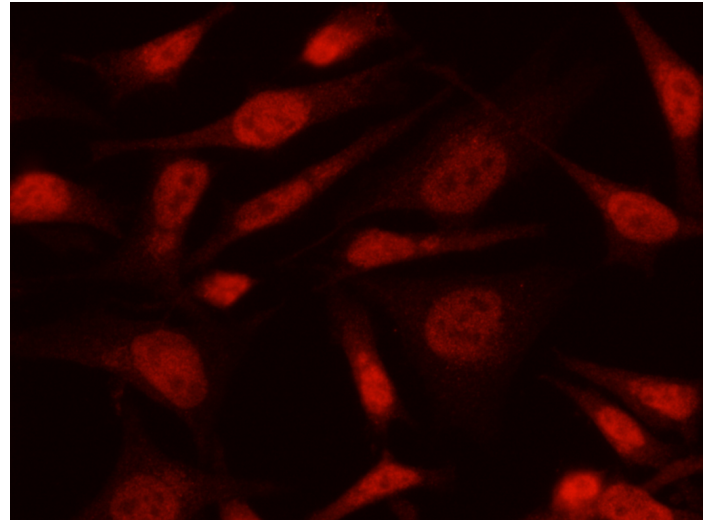
ENZ-ABS1036-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human MAX expression on Raji cells. The cells were treated according to manufacturer's manual (BD Pharmingen™ Cat. No. 554714), stained with purified anti-Human MAX, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of Human MAX in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-Human MAX monoclonal antibody (1:300) at 37° 1 hour. Then cells were stained with the Alexa Fluor® 594-conjugated goat Anti-rabbit IgG secondary antibody (red). Positive staining was localized to nucleus.

## Handling & Storage

**Handling** Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Protein max, MYC associated factor X

**Application** ELISA, Flow Cytometry, ICC, IF, IP, WB

**Clone** 103

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free. Carrier-free.

**Host** Rabbit

**Immunogen** Recombinant human MAX / MYC associated factor X.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation Dilutions/Conditions** ELISA (1:25'000-1:50'000)Flow Cytometry (1:100-1:500)Western Blot (1:500-1:1'000)Immunocytochemistry / Immunofluorescence (1:100-1:500)Immunoprecipitation (0.2-1 µL/mg of lysate)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**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-  
be@enzolifesciences.com](mailto:info-be@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)