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.

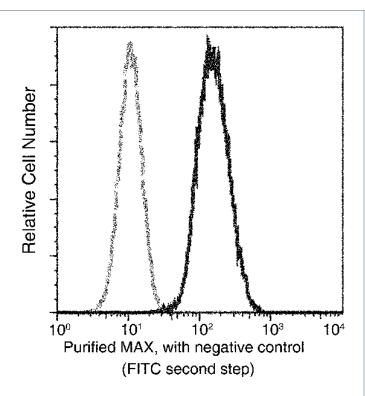
Ordering Information

Order Online »

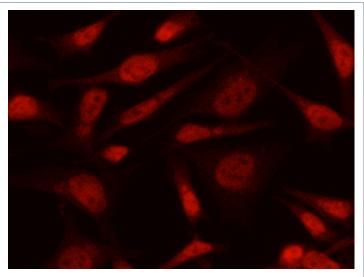
ENZ-ABS1036-0100 100μl

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, permeabilzed 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 μ L/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 infoLast modified: October 7, 2024

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

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com