

Bmf monoclonal antibody (12E10)

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-342-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Bcl-2-modifying factor

Application Flow Cytometry, IP, WB

Application Notes Excellent for Flow Cytometry.

Clone 12E10

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rat

Immunogen Recombinant mouse Bmf.

Isotype IgG1

Purity Detail Protein G-affinity purified.

Species Reactivity Human, Mouse

UniProt ID Q91ZE9

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



ENZO LIFE SCIENCES,
INC.
Phone: +1 914 361 6700
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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