

Bmf (mouse/rat) monoclonal antibody (17A9)

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-5080	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	ELISA, Flow Cytometry, IP, WB
Application Notes	Detects bands of ~25kDa and ~30kDa by Western blot, corresponding to endogenous Bmf and Bmf _L .
Clone	17A9
Crossreactivity	Does not cross-react with human Bmf.
Formulation	Liquid. Protein G affinity purified antibody in PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant full length mouse Bmf and synthetic peptides corresponding to aa 13-27 (D13VFQSEEDGEPGTQPG27) and aa 112-126 (A112GSPLGEPPEGQFL126) of mouse Bmf.
Isotype	IgG2a
Recommendation Dilutions/Conditions	Western Blot (1-2µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse, Rat
UniProt ID	Q91ZE9

Last modified: May 29, 2024



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)