

γ -Aminobutyric acid monoclonal antibody (5A9)

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

Ordering Information

[Order Online »](#)

BML-GA1160-0100	100 μ l
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GABA
Application	IHC, WB
Clone	5A9
Formulation	Liquid. In PBS solution containing 1.0% (w/v) BSA and 0.09% (w/v) sodium azide.
Host	Mouse
Immunogen	GABA conjugated to bovine serum albumin using glutaraldehyde.
Isotype	IgG1
Species Reactivity	Species independent
Specificity	Recognizes GABA. Species independent.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: February 14, 2025



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)