

Glial Fibrillary Acidic Protein monoclonal antibody (GA5)

GFAP is an intermediate filament protein expressed in astrocytes and other glial cells of the central nervous system. It serves as a widely used marker for astrocytic differentiation and CNS pathologies.

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

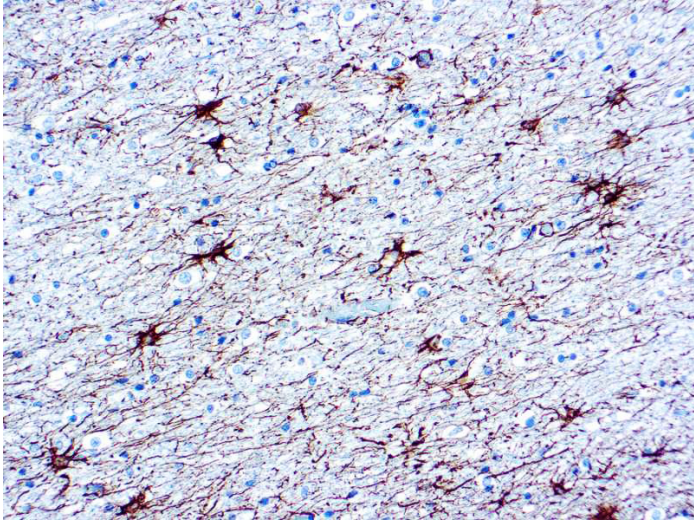
Ordering Information

[Order Online »](#)

ENZ-ABS1083-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human brain stained with GFAP antibody

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name GFAP

Application IHC (PS)

Application Notes Pretreatment – Citrate Buffer pH 6.0 (Manual/ Montage).
Incubation & Temperature 30 min @ RT

Clone GA5

Formulation Contains sodium azide as a preservative.

Host Mouse

Immunogen ALB/C mice were immunized with GFAP from porcine spinal cord.

Isotype IgG1κ

Purity Detail Purified immunoglobulin.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:50-1:200)

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](#).

Last modified: June 6, 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)