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

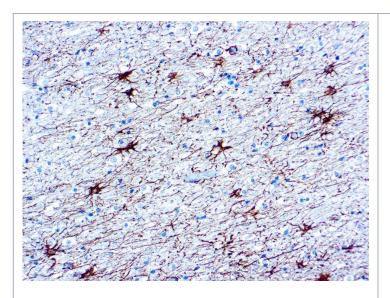
Order Online »

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.

IgG1ĸ

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



info-

eu@enzolifesciences.com