

GFAP Recombinant Monoclonal Mouse Antibody (rGA5) - Biotium Choice

A recombinant mouse monoclonal antibody that recognizes GFAP. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

GFAP Recombinant Monoclonal Mouse Antibody (rGA5) recognizes the glial fibrillary acidic protein (GFAP) expressed in neural tissues. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to a selection of CF® Dyes. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 7 bright and photostable CF® Dyes, including [near-IR CF®740](#)
- Suitable for immunohistochemistry, immunofluorescence, and western

GFAP is a 49-kDa type III intermediate filament protein found in neural tissues, serving as a marker that differentiates astrocytes from other glial cells during the development of the central nervous system. Although three splice variants of GFAP have been identified, the alpha isoform is the most abundantly expressed in astrocytes. GFAP is capable of co-assembling with vimentin and nestin in astrocytes, though these interactions are not essential for its filament formation. Similar to other intermediate filaments, GFAP assembly is regulated by phosphorylation and dephosphorylation of its N-terminal domain. Mutations in the GFAP gene have been linked to Alexander disease, and elevated GFAP expression has been observed in certain tumors derived from glial cells.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies – Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)

GFAP Recombinant Monoclonal Mouse Antibody (rGA5) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	1 mL	P046-488A-1ML	CF®488A Features
CF®568	562/584 nm	100 ug/mL	1 mL	P046-568-1ML	CF®568 Features
CF®594	593/615 nm	100 ug/mL	1 mL	P046-594-1ML	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	1 mL	P046-640R-1ML	CF®640R Features
CF®647	652/668 nm	100 ug/mL	1 mL	P046-647-1ML	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	1 mL	P046-680R-1ML	CF®680R Features
CF®740	742/767 nm	100 ug/mL	1 mL	P046-740-1ML	CF®740 Features

Product attributes

Antibody number	P046
Antibody reactivity (target)	GFAP
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary, Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rGA5
Isotype	IgG1
Molecular weight	50 kDa
Synonyms	GFAP; Glial fibrillary acidic protein
Human gene symbol	GFAP
SwissProt	P14136
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	Chicken, Human, Mouse, Pig, Rat
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user
Positive control	Brain, retina
Shipping condition	Room temperature
Storage Conditions	Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide
Antibody research areas	Neuroscience
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Expected antibody applications	IHC (FFPE) (published for clone), WB (published for clone)
Entrez gene ID	2670
Cell/tissue expression	Glia

This datasheet was generated on January 14, 2026 at 12:03:05 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/gfap-recombinant-monoclonal-mouse-antibody-rga5-biotium-choice/>