



Glowing products for science

## GFP Recombinant Monoclonal Mouse Antibody (r9F9.F9) – Biotium Choice

A recombinant mouse monoclonal antibody that recognizes the green fluorescent protein (GFP). The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

### Product Description

GFP (clone r9F9.F9) is a recombinant mouse monoclonal antibody that recognizes the green fluorescent protein (GFP). This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for fluorescence microscopy. The antibody is available purified, or conjugated to a selection of CF® Dyes or biotin.

### Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies Powered by CF® Dyes and Astral Leap™ Tandem Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes and Astral Leap™ tandem dyes, they are our top-recommended antibodies for flow cytometry and other applications.

#### Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for flow cytometry and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- Also available with [Astral Leap™ tandem dyes](#) for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

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



Call us : [800-304-5357](tel:800-304-5357)

#### Product attributes

Antibody reactivity (target)	GFP
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary, Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	r9F9.F9
Isotype	IgG1, kappa
Molecular weight	27 kDa
Synonyms	GFP, Green fluorescent protein
Verified antibody applications	IF (verified)
Antibody application notes	Immunofluorescence (HeLa): 1-5 ug/mL. Optimal concentration to be determined by end-user.
Positive control	HeLa or other human cell line
Shipping condition	Room temperature
Storage Conditions	Note: store unconjugated formats at -35 °C to -5 °C, Store at 2 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 research areas	Protein tag
Product origin	Animal
Filter by antibody application	IF (verified)
Antibody/conjugate formulation	Conjugates: PBS/50% glycerol/2 mg/mL rBSA/ 0.05% azide, Purified: PBS, no BSA, no azide

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF@405S</a>	411/431 nm	1 mg/mL	100 uL	<a href="#">P027-405S-100UL</a>	<a href="#">CF@405S Features</a>
<a href="#">CF@488A</a>	490/516 nm	1 mg/mL	100 uL	<a href="#">P027-488A-100UL</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	1 mg/mL	100 uL	<a href="#">P027-568-100UL</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@594</a>	593/615 nm	1 mg/mL	100 uL	<a href="#">P027-594-100UL</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@633</a>	629/650 nm	1 mg/mL	100 uL	<a href="#">P027-633-100UL</a>	<a href="#">CF@633 Features</a>
<a href="#">CF@640R</a>	642/663 nm	1 mg/mL	100 uL	<a href="#">P027-640R-100UL</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@647</a>	652/668 nm	1 mg/mL	100 uL	<a href="#">P027-647-100UL</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@660R</a>	662/682 nm	1 mg/mL	100 uL	<a href="#">P027-660R-100UL</a>	<a href="#">CF@660R Features</a>
<a href="#">CF@680</a>	681/698 nm	1 mg/mL	100 uL	<a href="#">P027-680-100UL</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@680R</a>	680/701 nm	1 mg/mL	100 uL	<a href="#">P027-680R-100UL</a>	<a href="#">CF@680R Features</a>
<a href="#">CF@740</a>	742/767 nm	1 mg/mL	100 uL	<a href="#">P027-740-100UL</a>	<a href="#">CF@740 Features</a>
<a href="#">CF@770</a>	770/797 nm	1 mg/mL	100 uL	<a href="#">P027-770-100UL</a>	<a href="#">CF@770 Features</a>
<a href="#">Biotin</a>	N/A	1 mg/mL	100 uL	<a href="#">P027-BTN-100UL</a>	
<a href="#">Purified (BSA-free)</a>	N/A	5 mg/mL	200 uL	<a href="#">P027-1MG</a>	
1000 uL	<a href="#">P027-5MG</a>				

This datasheet was generated on January 14, 2026 at 12:03:00 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/gfp-recombinant-monoclonal-mouse-antibody-r9f9-f9-biotium-choice/>