

# **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

	. 61	otlum	
00		-	
O.	Tax I	_0	
	Name March	-	_

Call us: 800-304-5357

Product attributes

Antibody reactivity (target)	GFP			
<b>Biotium Choice Antibody</b>	Primary			
Antibody type	Biotium Choice Flow Antibodies			
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 $^{\circ}$ C to -5 $^{\circ}$ C, Store at 2 to 8 $^{\circ}$ 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
CF®405S	411/431 nm	1 mg/mL	100 uL	P027-405S-100UL	CF®405S Features
CF®488A	490/516 nm	1 mg/mL	100 uL	P027-488A-100UL	CF®488A Features
CF®568	562/584 nm	1 mg/mL	100 uL	P027-568-100UL	CF®568 Features
CF®594	593/615 nm	1 mg/mL	100 uL	P027-594-100UL	CF®594 Features
CF®633	629/650 nm	1 mg/mL	100 uL	P027-633-100UL	CF®633 Features
CF®640R	642/663 nm	1 mg/mL	100 uL	P027-640R-100UL	CF®640R Features
CF®647	652/668 nm	1 mg/mL	100 uL	P027-647-100UL	CF®647 Features
CF®660R	662/682 nm	1 mg/mL	100 uL	P027-660R-100UL	CF®660R Features
CF®680	681/698 nm	1 mg/mL	100 uL	P027-680-100UL	CF®680 Features
CF®680R	680/701 nm	1 mg/mL	100 uL	P027-680R-100UL	CF®680R Features
CF®740	742/767 nm	1 mg/mL	100 uL	P027-740-100UL	CF®740 Features
CF®770	770/797 nm	1 mg/mL	100 uL	P027-770-100UL	CF®770 Features
<u>Biotin</u>	N/A	1 mg/mL	100 uL	P027-BTN-100UL	
Purified (BSA-free)	N/A	5 mg/mL	200 uL	P027-1MG	
1000 uL	P027-5MG				

This datasheet was generated on November 25, 2025 at 10:51:15 PM. 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/