

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](#)

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
Biotium Choice Antibody (SA-free)	488/508 nm	5 mg/mL	100 uL	P027-500B-000UL	CF®500B Features
			1000 uL	P027-5MG	

Product attributes

Antibody Number	P027
Antibody reactivity (target)	GFP
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice 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