

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.

[View our full catalog of Biotium Choice antibodies](#), featuring optimized and validated conjugates for microscopy or flow cytometry.

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

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
Biotium Choice (BSA-free)	519/599 nm	5 mg/mL	200 uL	P027-5MG-0001L	CF®500 Features
			1000 uL	P027-5MG	

This datasheet was generated on April 4, 2025 at 11:48:14 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/>