

## RFP Recombinant Monoclonal Mouse Antibody (r8E5.G7) - Biotium Choice



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

### Product Description

RFP Recombinant Monoclonal Mouse Antibody (r8E5.G7) is a recombinant mouse monoclonal antibody that recognizes the red fluorescent protein (RFP). 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 HRP.

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 bright and photostable CF® Dyes, including [near-IR CF®740](#), [CF®770](#), and [CF®790](#)
- Suitable for immunofluorescence and western

Red Fluorescent Protein (RFP) is isolated from sea anemone (*Discosoma sp.*, Uniprot ID: Q9U6Y8), and commonly used as a molecular tag to study protein localization, reporter gene expression, and protein-protein interactions. Although it exists as a homotetramer, monomeric variants have been constructed for use in various applications.

### 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](#)

## RFP Recombinant Monoclonal Mouse Antibody (r8E5.G7) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF@488A</a>	490/516 nm	100 ug/mL	500 uL	<a href="#">P053-488A-500UL</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	100 ug/mL	500 uL	<a href="#">P053-568-500UL</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@594</a>	593/615 nm	100 ug/mL	500 uL	<a href="#">P053-594-500UL</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@640R</a>	642/663 nm	100 ug/mL	500 uL	<a href="#">P053-640R-500UL</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@647</a>	652/668 nm	100 ug/mL	500 uL	<a href="#">P053-647-500UL</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@680</a>	680/701 nm	100 ug/mL	500 uL	<a href="#">P053-680-500UL</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@680R</a>	680/701 nm	100 ug/mL	500 uL	<a href="#">P053-680R-500UL</a>	<a href="#">CF@680R Features</a>
<a href="#">CF@740</a>	742/767 nm	100 ug/mL	500 uL	<a href="#">P053-740-500UL</a>	<a href="#">CF@740 Features</a>
<a href="#">CF@770</a>	770/797 nm	100 ug/mL	500 uL	<a href="#">P053-770-500UL</a>	<a href="#">CF@770 Features</a>
<a href="#">CF@790</a>	783/808 nm	100 ug/mL	500 uL	<a href="#">P053-790-500UL</a>	<a href="#">CF@790 Features</a>
<a href="#">HRP</a>		100 ug/mL	500 uL	<a href="#">P053-HRP-500UL</a>	

### Product attributes

Expected antibody applications	WB (published)
Antibody number	P053
Antibody reactivity (target)	RFP
Synonyms	sRed2; mRFP; red fluorescent; Red fluorescent protein; RFP; RFP tag
<b>Biotium Choice Antibody</b>	Primary, Tag Antibody
Antibody type	Biotium Choice, Primary, Tag Antibody
Host species	Mouse
Clonality	Monoclonal, Recombinant Monoclonal
Clone	r8E5.G7
Isotype	IgG2a
Verified antibody applications	IF (verified)
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user
Shipping condition	Room temperature
Storage Conditions	Store CF® Dye conjugates at 2 °C to 8 °C. Store HRP conjugates at -10 °C to -35 °C, Protect 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	CF® Dye Conjugates: PBS/2 mg/mL rBSA/ 0.05% sodium azide, HRP conjugates: PBS/50% glycerol/2 mg/mL rBSA
Antibody research areas	Protein tag
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line

This datasheet was generated on June 3, 2026 at 05:55:58 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rfp-recombinant-monoclonal-mouse-antibody-r8e5-g7-biotium-choice/>