



Glowing products for science

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

Fluorescein (FITC) Recombinant Monoclonal Mouse Antibody (r1F8.1E4) - Biotium Choice



Recombinant Monoclonal mouse anti-fluorescein (FITC) antibody suitable for IF, flow, and Western blot. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

Product Description

Recombinant Monoclonal mouse anti-fluorescein (FITC) antibody labeled with our superior CF® dyes.

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](#) and [CF®790](#)
- Suitable for immunofluorescence, IHC, western, and flow cytometry

FITC (fluorescein isothiocyanate) is a fluorochrome dye that absorbs ultraviolet or blue light causing molecules to become excited and emit a visible yellow-green light. This emission ceases upon removal of the light causing the excitation. Fluorochrome labeling provides rapid, accurate localization of antigen-antibody interaction when one of the reactants is part of a cell, tissue or other biological structure. FITC is a commonly used marker for antibodies in immunofluorescent techniques since the conjugation of FITC to proteins is relatively easy and does not, in general, destroy the biological activity of the labeled protein. FITC is widely used as a hapten to label different proteins. Anti-FITC antibodies can be used for isolating FITC labeled cells by magnetic microbeads based sorting, as well as removal of background signal.

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](#)

Fluorescein (FITC) Recombinant Monoclonal Mouse Antibody (r1F8.1E4) - Biotium Choice

Conjugation	Ex/Em	Size	Catalog No.
CF®405M	408/452 nm	500 uL	P063-405M-500UL
CF®488A	490/515 nm	500 uL	P063-488A-500UL
CF®568	562/584 nm	500 uL	P063-568-500UL
CF®594	593/614 nm	500 uL	P063-594-500UL
CF®640R	642/662 nm	500 uL	P063-640R-500UL
CF®647	652/668 nm	500 uL	P063-647-500UL
CF®680R	680/701 nm	500 uL	P063-680R-500UL
CF®740	742/767 nm	500 uL	P063-740-500UL
CF®790	783/808 nm	500 uL	P063-790-500UL

This datasheet was generated on September 22, 2026 at 02:33:42 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/fluorescein-fitc-recombinant-monoclonal-mouse-antibody-r1f8-1e4-biotium-choice/>

Product attributes

Antibody number	P063
Antibody reactivity (target)	Fluorescein (FITC)
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice, Primary
Host species	Mouse
Clonality	Monoclonal, Recombinant Monoclonal
Clone	r1F8.1E4
Isotype	IgG2, kappa
Synonyms	Fluorescein; FITC
Verified antibody applications	IF (verified)
Expected antibody applications	Flow (intracellular), Flow (surface), IF, WB
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, 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
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Volume per assay	5 uL/test