



Glowing products for science

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

## Fluorescein (FITC) Monoclonal Mouse Antibody (1F8.1E4)

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



### Product Description

This is a monoclonal mouse anti-fluorescein (FITC) antibody labeled with our superior CF® dyes.

- Available in 6 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

See our full selection of [anti-tag and anti-hapten antibody conjugates](#).

**Note:** Conjugates of blue fluorescent dyes like CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

### Monoclonal Mouse Anti-Fluorescein (FITC)

| Conjugation      | Ex/Em                        | Size           | Catalog No.                |
|------------------|------------------------------|----------------|----------------------------|
| CF®405M          | 408/452 nm                   | 50 uL (100 ug) | <a href="#">20214-1</a>    |
| 0.25 mL (500 ug) | <a href="#">20214</a>        |                |                            |
| CF®488A          | 490/515 nm                   | 50 uL (100 ug) | <a href="#">20210-1</a>    |
| 0.25 mL (500 ug) | <a href="#">20210</a>        |                |                            |
| CF®594           | 593/614 nm                   | 50 uL (100 ug) | <a href="#">20211-1</a>    |
| 0.25 mL (500 ug) | <a href="#">20211</a>        |                |                            |
| CF®633           | 630/650 nm                   | 50 uL (100 ug) | <a href="#">20212-1</a>    |
| 0.25 mL (500 ug) | <a href="#">20212</a>        |                |                            |
| CF®640R          | 642/662 nm                   | 50 uL (100 ug) | <a href="#">20213-1</a>    |
| 0.25 mL (500 ug) | <a href="#">20213</a>        |                |                            |
| CF®660R          | 663/682 nm                   | 50 uL (100 ug) | <a href="#">20399-50uL</a> |
| 0.25 mL (500 ug) | <a href="#">20399-0.25mL</a> |                |                            |

View our full selection of [Secondary Antibodies](#), or search our catalog using our [Antibody Finder](#). Alternatively, you can view our [secondary antibody product listings](#) with catalog numbers.

CF® Dyes offer exceptional brightness and photostability. For more information see our [CF® Dye technology page](#).

#### Storage and Handling

**Liquid format:** Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Liquid format antibodies contain 50% glycerol and will not freeze at -20°C.

**Lyophilized format:** Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Reconstitute antibodies in water using the indicated volumes below:

CF® Dye and biotin conjugates: add 0.5 mL dH<sub>2</sub>O

HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O

Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store reconstituted antibody at -20°C and protect from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20°C.

Optional: A preservative such as 0.05% sodium azide (final concentration) can be added to CF® Dye and biotin conjugates. Do not add sodium azide to HRP conjugates.

**Note:** Storage of the antibody for more than a day at final working dilution is not recommended.

CF is a registered trademark of Biotium, Inc.

### References

Download a list of [CF® dye references](#).

#### Product attributes

|                                     |   |
|-------------------------------------|---|
| Antibody type                       | Tag Antibody  |
| Antibody reactivity (target)        | FITC  |
| Host species                        | Mouse   |
| Clonality                           | Monoclonal  |
| Clone                               | 1F8.1E4   |
| Isotype                             | IgG1, kappa   |
| Antibody/conjugate formulation      | Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide  |
| Concentration                       | 2 mg/mL   |
| Secondary/tag antibody applications | Flow cytometry, IHC, IF (cells or tissue sections), Western blot  |
| Product origin                      | Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |

This datasheet was generated on December 15, 2025 at 07:46:08 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-mouse-anti-fitc-antibody/>