

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. 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 dH₂O HRP or DNP conjugates: add 1 mL dH₂O 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.

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

References

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

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