



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: techsupport@biotium.com

GFP Monoclonal Mouse Antibody (9F9.F9)

Monoclonal mouse anti-GFP antibody labeled with our superior CF® dyes.



Product Description

This is a monoclonal mouse anti-GFP (green fluorescent protein) antibody labeled with our superior CF® Dyes.

- Available in 9 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Biotium now offers a [Recombinant GFP Monoclonal Mouse \(r9F9.F9\) Biotium Choice Antibody](#) which offers comparable performance at a lower cost.

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

Product attributes

Antibody type	Tag Antibody
Antibody reactivity (target)	GFP
Host species	Mouse
Clonality	Monoclonal
Clone	9F9.F9
Isotype	IgG1, kappa
Immunogen	Fusion protein corresponding to Aequorea victoria full length GFP, aa 1-246
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Concentration	1 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.

GFP Monoclonal Mouse Antibody

Conjugation	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	100 uL (100 ug)	20215
CF®568	562/583 nm	100 uL (100 ug)	20480
CF®594	593/614 nm	100 uL (100 ug)	20216
CF®633	630/650 nm	100 uL (100 ug)	20217
CF®640R	642/662 nm	100 uL (100 ug)	20218
CF®660R	663/682 nm	100 uL (100 ug)	20481
CF®680	681/698 nm	100 uL (100 ug)	20219
CF®680R	680/701 nm	100 uL (100 ug)	20482
CF®770	770/797 nm	100 uL (100 ug)	20220

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.

References

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

This datasheet was generated on September 9, 2025 at 06:04:34 PM. Visit product page to check for updated information before use.

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