



Glowing products for science

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

## Goat Anti-Mouse IgM (Mu Chain)

Goat anti-mouse IgM ( $\mu$  chain) secondary antibody labeled with our superior CF® dyes.



### Product Description

This is a goat anti-mouse IgM ( $\mu$  chain) secondary antibody labeled with our bright and photostable CF® dyes. The antibodies react specifically with IgM heavy chains ( $\mu$  chains) and not with immunoglobulin light chains.

- Available in 5 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology

#### Goat Anti-Mouse IgM (Mu Chain), CF® Dye Conjugates

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@488A</a>	490/515 nm	50 $\mu$ L (100 $\mu$ g)	<a href="#">20840-50uL</a>	<a href="#">CF@488A Features</a>
		0.25mL (500 $\mu$ g)	<a href="#">20840-250uL</a>	
<a href="#">CF@555</a>	555/565 nm	50 $\mu$ L (100 $\mu$ g)	<a href="#">20485-50uL</a>	<a href="#">CF@555 Features</a>
		0.25mL (500 $\mu$ g)	<a href="#">20485-250uL</a>	
<a href="#">CF@680</a>	681/698 nm	50 $\mu$ L (100 $\mu$ g)	<a href="#">20384-50uL</a>	<a href="#">CF@680 Features</a>
		0.25mL (500 $\mu$ g)	<a href="#">20384-250uL</a>	
<a href="#">CF@680R</a>	680/701 nm	50 $\mu$ L (100 $\mu$ g)	<a href="#">20841-50uL</a>	<a href="#">CF@680R Features</a>
		0.25mL (500 $\mu$ g)	<a href="#">20841-250uL</a>	
<a href="#">CF@770</a>	770/797 nm	50 $\mu$ L (100 $\mu$ g)	<a href="#">20385-50uL</a>	<a href="#">CF@770 Features</a>
		0.25mL (500 $\mu$ g)	<a href="#">20385-250uL</a>	

View our full selection of bright and specific [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 curated [CF® Dye references](#).

This datasheet was generated on August 29, 2025 at 06:19:39 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-mouse-igm-mu-chain-2-mgml/>

#### Product attributes

Clonality	Polyclonal
Antibody type	Secondary
Concentration	2 mg/mL
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
Species reactivity	Mouse
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.
Host species	Goat
Antibody reactivity (target)	Mouse IgM