

## Goat Anti-Mouse IgM (Mu Chain)

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



### Product Description

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

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

### 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 ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Host species	Goat
Antibody reactivity (target)	Mouse IgM

### Goat Anti-Mouse IgM (Mu Chain), CF<sup>®</sup> 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<sup>®</sup> Dyes offer exceptional brightness and photostability. For more information see our [CF<sup>®</sup> 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<sup>®</sup> 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<sup>®</sup> 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<sup>®</sup> Dye references](#).

This datasheet was generated on May 18, 2025 at 03:26:48 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/>