

## Goat Anti-Mouse IgM (Mu Chain)

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



### 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  |

## 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><br>0.25mL (500 ug) | 490/515 nm<br><a href="#">20840-250uL</a> | 50 uL (100 ug) | <a href="#">20840-50uL</a> | <a href="#">CF@488A Features</a> |
| <a href="#">CF@555</a><br>0.25mL (500 ug)  | 555/565 nm<br><a href="#">20485-250uL</a> | 50 uL (100 ug) | <a href="#">20485-50uL</a> | <a href="#">CF@555 Features</a>  |
| <a href="#">CF@680</a><br>0.25mL (500 ug)  | 681/698 nm<br><a href="#">20384-250uL</a> | 50 uL (100 ug) | <a href="#">20384-50uL</a> | <a href="#">CF@680 Features</a>  |
| <a href="#">CF@680R</a><br>0.25mL (500 ug) | 680/701 nm<br><a href="#">20841-250uL</a> | 50 uL (100 ug) | <a href="#">20841-50uL</a> | <a href="#">CF@680R Features</a> |
| <a href="#">CF@770</a><br>0.25mL (500 ug)  | 770/797 nm<br><a href="#">20385-250uL</a> | 50 uL (100 ug) | <a href="#">20385-50uL</a> | <a href="#">CF@770 Features</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 June 3, 2026 at 03:28:53 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/>