

Biotin Monoclonal Mouse Antibody (3D6.6)

Monoclonal mouse anti-biotin antibody labeled with our superior CF® Dyes.

Blottum

Product Description

This is a monoclonal mouse anti-biotin antibody labeled with our superior CF® Dyes. The antibody is useful for detecting biotin conjugated to antibodies or other proteins.

- Available in 8 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®405S 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.

Product attributes

Call us: 800-304-5357

| Antibody type | Tag Antibody |
|-------------------------------------|---|
| Antibody reactivity (target) | Biotin |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | 3D6.6 |
| Isotype | IgG1, kappa |
| Concentration | 2 mg/mL |
| Antibody/conjugate formulation | Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide |
| Storage Conditions | Store at -10 to -35 °C, Protect from light |
| 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. |
| | |

Email: techsupport@biotium.com

Monoclonal Mouse Anti-Biotin

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|-------------|------------|------------------|----------------|------------------|
| CF®405S | 404/431 nm | 50 uL (100 ug) | <u>20203-1</u> | CF®405S Features |
| | | 0.25 mL (500 ug) | <u>20203</u> | |
| CF®488A | 490/515 nm | 50 uL (100 ug) | <u>20204-1</u> | CF®488A Features |
| | | 0.25 mL (500 ug) | <u>20204</u> | |
| CF®568 | 562/584 nm | 50 uL (100 ug) | <u>20502-1</u> | CF®568 Features |
| | | 0.25 mL (500 ug) | <u>20502</u> | |
| CF®594 | 593/614 nm | 50 uL (100 ug) | <u>20205-1</u> | CF®594 Features |
| | | 0.25 mL (500 ug) | <u>20205</u> | |
| CF®633 | 630/650 nm | 50 uL (100 ug) | <u>20206-1</u> | CF®633 Features |
| | | 0.25 mL (500 ug) | <u>20206</u> | |
| CF®640R | 642/662 nm | 50 uL (100 ug) | <u>20207-1</u> | CF®640R Features |
| | | 0.25 mL (500 ug) | <u>20207</u> | |
| CF®750 | 755/777 nm | 50 uL (100 ug) | 20501-50uL | CF®750 Features |
| | | 0.25 mL (500 ug) | 20501-250uL | |
| CF®770 | 770/797 nm | 50 uL (100 ug) | 20367-50uL | CF®770 Features |
| | | 0.25 mL (500 ug) | 20367-250uL | |

View our full selection of <u>Secondary Antibodies</u>, or search our catalog using our <u>Antibody Finder</u>. Alternatively, you can view our <u>secondary antibody product listings</u> 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 CF® dye references.

This datasheet was generated on October 30, 2025 at 12:15:36 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-mouse-anti-biotin-antibody/