

## Biotin Monoclonal Mouse Antibody (3D6.6)

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



### Product attributes

<b>Antibody type</b>	Primary, Tag Antibody
<b>Antibody reactivity (target)</b>	Biotin
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	3D6.6
<b>Isotype</b>	IgG1, kappa
<b>Concentration</b>	2 mg/mL (CF® dye conjugates), 0.1 mg/mL (HRP conjugate)
<b>Antibody/conjugate formulation</b>	HRP conjugates: PBS/50% glycerol/2 mg/mL rBSA, Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
<b>Storage Conditions</b>	Store at -10 to -35 °C, Protect from light
<b>Secondary/tag antibody applications</b>	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
<b>Product origin</b>	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.

### 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, and HRP
- 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.

### Monoclonal Mouse Anti-Biotin

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®405S	404/431 nm	50 uL (100 ug)	<a href="#">20203-1</a>	<a href="#">CF®405S Features</a>
250 uL (500 ug)	<a href="#">20203</a>			
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20204-1</a>	<a href="#">CF®488A Features</a>
250 uL (500 ug)	<a href="#">20204</a>			
CF®568	562/584 nm	50 uL (100 ug)	<a href="#">20502-1</a>	<a href="#">CF®568 Features</a>
250 uL (500 ug)	<a href="#">20502</a>			
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20205-1</a>	<a href="#">CF®594 Features</a>
250 uL (500 ug)	<a href="#">20205</a>			
CF®633	630/650 nm	50 uL (100 ug)	<a href="#">20206-1</a>	<a href="#">CF®633 Features</a>
250 uL (500 ug)	<a href="#">20206</a>			
CF®640R	642/662 nm	50 uL (100 ug)	<a href="#">20207-1</a>	<a href="#">CF®640R Features</a>
250 uL (500 ug)	<a href="#">20207</a>			
CF®750	755/777 nm	50 uL (100 ug)	<a href="#">20501-50UL</a>	<a href="#">CF®750 Features</a>
250 uL (500 ug)	<a href="#">20501-250UL</a>			
CF®770	770/797 nm	50 uL (100 ug)	<a href="#">20367-50UL</a>	<a href="#">CF®770 Features</a>
250 uL (500 ug)	<a href="#">20367-250UL</a>			
HRP	N/A	500 uL (1 mg)	<a href="#">20503-500UL</a>	

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 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 [CF® dye references](#).

This datasheet was generated on June 3, 2026 at 04:38:15 PM. Visit product page to check for updated information before use.

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