

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

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



### 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. 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.

### Product attributes

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

### References

Download a list of [CF® dye references](#).

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>
		0.25 mL (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>
		0.25 mL (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>
		0.25 mL (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>
		0.25 mL (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>
		0.25 mL (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>
		0.25 mL (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>
		0.25 mL (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>
		0.25 mL (500 ug)	<a href="#">20367-250uL</a>	