

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.

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
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.

Monoclonal Mouse Anti-Biotin

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®405S	404/431 nm	50 uL (100 ug)	20203-1	CF®405S Features
		0.25 mL (500 ug)	20203	
CF®488A	490/515 nm	50 uL (100 ug)	20204-1	CF®488A Features
		0.25 mL (500 ug)	20204	
CF®568	562/584 nm	50 uL (100 ug)	20502-1	CF®568 Features
		0.25 mL (500 ug)	20502	
CF®594	593/614 nm	50 uL (100 ug)	20205-1	CF®594 Features
		0.25 mL (500 ug)	20205	
CF®633	630/650 nm	50 uL (100 ug)	20206-1	CF®633 Features
		0.25 mL (500 ug)	20206	
CF®640R	642/662 nm	50 uL (100 ug)	20207-1	CF®640R Features
		0.25 mL (500 ug)	20207	
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 [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 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 September 2, 2025 at 01:07:15 AM. Visit product page to check for updated information before use.

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