

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 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.

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) 0.25 mL (500 ug)	20203-1 20203	CF®405S Features
CF®488A	490/515 nm	50 uL (100 ug) 0.25 mL (500 ug)	20204-1 20204	CF®488A Features
CF®568	562/584 nm	50 uL (100 ug) 0.25 mL (500 ug)	20502-1 20502	CF®568 Features
CF®594	593/614 nm	50 uL (100 ug) 0.25 mL (500 ug)	20205-1 20205	CF®594 Features
CF®633	630/650 nm	50 uL (100 ug) 0.25 mL (500 ug)	20206-1 20206	CF®633 Features
CF®640R	642/662 nm	50 uL (100 ug) 0.25 mL (500 ug)	20207-1 20207	CF®640R Features
CF®750	755/777 nm	50 uL (100 ug) 0.25 mL (500 ug)	20501-50uL 20501-250uL	CF®750 Features
CF®770	770/797 nm	50 uL (100 ug) 0.25 mL (500 ug)	20367-50uL 20367-250uL	CF®770 Features