

6X His Tag Monoclonal Mouse Antibody

Monoclonal mouse anti-6X His Tag antibody labeled with our superior CF® dyes.



Product Description

This is a monoclonal mouse anti-6X His Tag antibody labeled with our superior CF® dyes. The antibody recognizes the 6X His epitope tag (His-His-His-His-His-His) fused to either the amino- or carboxy- termini of targeted proteins.

- Available in 5 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](#). 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)	6X His Tag
Host species	Mouse
Clonality	Monoclonal
Clone	33D10.D2.G8
Isotype	IgG1, kappa
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Concentration	1 mg/mL
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.
CF®488A	490/515 nm	50 uL (50 ug)	20228
CF®594	593/614 nm	50 uL (50 ug)	20229
CF®640R	642/662 nm	50 uL (50 ug)	20237
CF®680	681/698 nm	50 uL (50 ug)	20359
CF®770	770/797 nm	50 uL (50 ug)	20360