

## 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](#).

### Product attributes

<b>Antibody type</b>	Primary, Tag Antibody
<b>Antibody reactivity (target)</b>	6X His Tag
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	33D10.D2.G8
<b>Isotype</b>	IgG1, kappa
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
<b>Concentration</b>	1 mg/mL
<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.

## Monoclonal Mouse Anti-6X His

Conjugation	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	50 uL (50 ug)	<a href="#">20228</a>
CF®594	593/614 nm	50 uL (50 ug)	<a href="#">20229</a>
CF®640R	642/662 nm	50 uL (50 ug)	<a href="#">20237</a>
CF®680	681/698 nm	50 uL (50 ug)	<a href="#">20359</a>
CF®770	770/797 nm	50 uL (50 ug)	<a href="#">20360</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 04:05:35 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-mouse-anti-6x-his-antibody/>