

## 6X His Tag Recombinant Monoclonal Mouse Antibody (2405.H6) - Biotium Choice



Recombinant monoclonal mouse anti-6X His Tag antibody labeled with our superior CF® Dyes.

### Product Description

6X His Tag (clone 2405.H6) is a recombinant monoclonal mouse antibody that recognizes the 6X His epitope tag (His-His-His-His-His-His) fused to either the amino- or carboxy- termini of targeted proteins. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to a selection of CF® Dyes and other conjugates.

- Available in bright and photostable CF® Dyes, as well as RPE, APC, and HRP
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

See our full selection of [anti-tag and anti-hapten antibody conjugates](#).

### Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies Powered by CF® Dyes and Astral Leap™ Tandem Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes and Astral Leap™ tandem dyes, they are our top-recommended antibodies for flow cytometry and other applications.

#### Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for flow cytometry and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- Also available with [Astral Leap™ tandem dyes](#) for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)

## 6X His Tag Recombinant Monoclonal Mouse Antibody (2405.H6) – Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF®405S</a>	411/431 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-405S-100UL</a>	<a href="#">CF®405S Features</a>
<a href="#">CF®488A</a>	490/516 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-488A-100UL</a>	<a href="#">CF®488A Features</a>
<a href="#">CF®568</a>	562/584 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-568-100UL</a>	<a href="#">CF®568 Features</a>
<a href="#">CF®594</a>	593/615 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-594-100UL</a>	<a href="#">CF®594 Features</a>
<a href="#">CF®640R</a>	642/663 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-640R-100UL</a>	<a href="#">CF®640R Features</a>
<a href="#">CF®647</a>	652/668 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-647-100UL</a>	<a href="#">CF®647 Features</a>
<a href="#">CF®680</a>	681/698 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-680-100UL</a>	<a href="#">CF®680 Features</a>
<a href="#">CF®680R</a>	680/701 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-680R-100UL</a>	<a href="#">CF®680R Features</a>
<a href="#">CF®740</a>	742/767 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-740-100UL</a>	<a href="#">CF®740 Features</a>
<a href="#">CF®770</a>	770/797 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-770-100UL</a>	<a href="#">CF®770 Features</a>
<a href="#">CF®790</a>	783/808 nm	1 mg/mL	100 uL (100 ug)	<a href="#">P024-790-100UL</a>	<a href="#">CF®790 Features</a>
<a href="#">RPE</a>	N/A	0.5 mg/mL	100 uL (50 ug)	<a href="#">P024-RPE-100UL</a>	
<a href="#">APC</a>	N/A	0.5 mg/mL	100 uL (50 ug)	<a href="#">P024-APC-100UL</a>	
<a href="#">HRP</a>	N/A	0.5 mg/mL	100 uL (50 ug)	<a href="#">P024-HRP-100UL</a>	

### Product attributes

Antibody number	#P024
Biotium Choice Antibody	Primary, Tag Antibody
Antibody type	Biotium Choice, Primary, Tag Antibody
Clonality	Monoclonal, Recombinant Monoclonal
Host species	Mouse
Clone	2405.H6
Isotype	IgG1
Antibody reactivity (target)	6X His Tag
Synonyms	Histidine tag, 6 His epitope tag, 6-His Tag, 6X His, 6x-His Tag, 6X-His, H, Hexa His tag, HHHHHH epitope tag, HH-HHHH tag, His Tag, HIS; His6-Tag, polyhistidine tag, polyHistidine, Poly-histidine
Species reactivity	Bovine, Chicken, Duck, Human, Monkey, Mouse, Pig, Rat, Yeast
Verified antibody applications	IF (verified), WB (verified), Near-IR WB (verified)
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Positive control	HeLa or other human cell line
Antibody application notes	Immunofluorescence (HeLa): 1-5 ug/mL, Near-IR WB: 0.1-0.5 ug/mL. Optimal concentration to be determined by end-user.
Antibody research areas	Protein tag
Antibody/conjugate formulation	Conjugates: PBS/50% glycerol/2 mg/mL rBSA/ 0.05% azide, HRP conjugates: PBS/50% glycerol/2 mg/mL rBSA
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light
Product origin	Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line; recombinant BSA produced in Chinese hamster ovary cells.
Regulatory status	For research use only (RUO)
Shipping condition	Room temperature
Expected antibody applications	IF
Secondary/tag antibody applications	IF (cells or tissue sections), Western blot

### References

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

This datasheet was generated on June 3, 2026 at 04:05:31 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/6x-his-tag-recombinant-monoclonal-mouse-antibody-2405-h6-biotium-choice/>