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 10 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 <u>Astral Leap™ tandem dyes</u> for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

View our full catalog of Biotium Choice antibodies

Product attributes

applications

Antibody number	#P024			
Biotium Choice Antibody	Tag Antibody			
Antibody type	Biotium Choice Tag Antibodies, Tag Antibody			
Clonality	Recombinant Monoclonal			
Host species	Mouse			
Clone	2405.H6			
Isotype	lgG1			
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, HHHHHHH epitope tag, HHHHHHH 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)			
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.			
D	For research use only (RUO)			
Regulatory status	roi researcii use only (noo)			
Shipping condition	Room temperature			
• ,	,, ,			

Call us: 800-304-5357 Email: btinfo@biotium.com

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

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

References

Download a list of CF® Dye references.

This datasheet was generated on November 2, 2025 at 03:15:37 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/