

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.

- Available in 10 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](#). Also view our line of highly optimized [Biotium Choice Antibodies](#) for flow cytometry.

CF® Dyes offer exceptional brightness and photostability. For more information see our [CF® Dye technology page](#).

Product attributes

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	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)
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

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
CF@488A	490/516 nm	1 mg/mL	100 uL (100 ug)	P024-488A-100UL	CF@488A Features
CF@568	562/584 nm	1 mg/mL	100 uL (100 ug)	P024-568-100UL	CF@568 Features
CF@594	593/615 nm	1 mg/mL	100 uL (100 ug)	P024-594-100UL	CF@594 Features
CF@640R	642/663 nm	1 mg/mL	100 uL (100 ug)	P024-640R-100UL	CF@640R Features
CF@647	652/668 nm	1 mg/mL	100 uL (100 ug)	P024-647-100UL	CF@647 Features
CF@680	681/698 nm	1 mg/mL	100 uL (100 ug)	P024-680-100UL	CF@680 Features
CF@680R	680/701 nm	1 mg/mL	100 uL (100 ug)	P024-680R-100UL	CF@680R Features
CF@740	742/767 nm	1 mg/mL	100 uL (100 ug)	P024-740-100UL	CF@740 Features
CF@770	770/797 nm	1 mg/mL	100 uL (100 ug)	P024-770-100UL	CF@770 Features

References

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

This datasheet was generated on January 15, 2025 at 05:49:49 AM. 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/>