



Glowing products for science

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 [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. | |
|-------------------------|------------|-----------|-----------------|---------------------------------|----------------------------------|
| 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 |
| RPE | N/A | 0.5 mg/mL | 100 uL (50 ug) | P024-RPE-100UL | |
| APC | 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 December 22, 2025 at 11:23:11 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/>



Call us : [800-304-5357](tel:800-304-5357)

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, HHHHHH 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 |