

HA Tag Recombinant Monoclonal Mouse Antibody (r12CA5) - Biotium Choice



Recombinant monoclonal mouse anti-HA Tag antibody labeled with our superior CF® Dyes. A recombinant mouse monoclonal antibody that recognizes the HA (hemagglutinin) epitope tag. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

Product Description

HA Tag (clone r12CA5) is a recombinant monoclonal mouse antibody that recognizes the HA (hemagglutinin) epitope tag. 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 HRP.

- Available in bright and photostable CF® Dyes and HRP
- Suitable for western and immunofluorescence

The HA tag (hemagglutinin tag) is an epitope tag widely used to localize gene products across diverse cell types, investigate the topology of proteins and protein complexes, identify associated proteins, and characterize newly discovered, low-abundance, or poorly immunogenic proteins in situations where protein-specific antibodies are unavailable. Derived from the influenza surface glycoprotein hemagglutinin, the HA tag is routinely co-expressed with proteins of interest to facilitate their identification and purification for downstream functional analysis.

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](#)

HA Tag Recombinant Monoclonal Mouse Antibody (r12CA5) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF@488A	490/516 nm	100 ug/mL	500 uL (50 ug)	P054-488A-500UL	CF@488A Features
CF@568	562/584 nm	100 ug/mL	500 uL (50 ug)	P054-568-500UL	CF@568 Features
CF@594	593/615 nm	100 ug/mL	500 uL (50 ug)	P054-594-500UL	CF@594 Features
CF@640R	642/663 nm	100 ug/mL	500 uL (50 ug)	P054-640R-500UL	CF@640R Features
CF@647	652/668 nm	100 ug/mL	500 uL (50 ug)	P054-647-500UL	CF@647 Features
CF@680	681/698 nm	100 ug/mL	500 uL (50 ug)	P054-680-500UL	CF@680 Features
CF@680R	680/701 nm	100 ug/mL	500 uL (50 ug)	P054-680R-500UL	CF@680R Features
CF@740	742/767 nm	100 ug/mL	500 uL (50 ug)	P054-740-500UL	CF@740 Features
CF@770	770/797 nm	100 ug/mL	500 uL (50 ug)	P054-770-500UL	CF@770 Features
CF@790	783/808 nm	100 ug/mL	500 uL (50 ug)	P054-790-500UL	CF@790 Features
HRP	N/A	100 ug/mL	500 uL (50 ug)	P054-HRP-500UL	

Product attributes

Antibody number	#P054
Biotium Choice Antibody	Primary, Tag Antibody
Antibody type	Biotium Choice, Primary, Tag Antibody
Clonality	Monoclonal, Recombinant Monoclonal
Host species	Mouse
Clone	r12CA5
Isotype	IgG2b
Antibody reactivity (target)	HA Tag
Synonyms	HA epitope tag, hemagglutinin, YPYDVPDYA, Hemagglutinin tag
Immunogen	9 amino acid peptide to influenza virus hemagglutinin
Species reactivity	Bovine, Chicken, Duck, Human, Monkey, Mouse, Pig, Rat, Yeast
Verified antibody applications	IF (verified), WB (verified)
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody application notes	Immunofluorescence: 0.25-1 ug/mL, Optimal concentration to be determined by end-user, Western Blot (fluorescent): 100-500 ng/mL, Western Blot (HRP): 10-50 ng/mL
Antibody research areas	Protein tag
Antibody/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide, HRP conjugates: PBS/50% glycerol/2 mg/mL rBSA
Storage Conditions	Store CF® Dye conjugates at 2 °C to 8 °C, Store HRP conjugates at -10 °C to -35 °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	IHC (FFPE) (published for clone), IP (published for clone)
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 04:53:26 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ha-tag-recombinant-monoclonal-mouse-antibody-r12ca5-biotium-choice/>