

c-Fos Recombinant Monoclonal Mouse Antibody (r2H2) - Biotium Choice

A recombinant mouse monoclonal antibody that recognizes FBJ murine osteosarcoma viral oncogene homolog, commonly known as Cellular protooncogene Fos (c-Fos). This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

c-Fos Recombinant Monoclonal Mouse Antibody (r2H2) is a recombinant mouse monoclonal antibody that recognizes FBJ murine osteosarcoma viral oncogene homolog, commonly known as Cellular protooncogene Fos (c-Fos). 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. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 7 bright and photostable CF® Dyes, including [near-IR CF®740](#)
- Suitable for immunohistochemistry, immunofluorescence, and western

c-Fos, encoded by the C-FOS gene on chromosome 14, is a member of the Fos transcription factor family and forms the AP-1 complex with c-Jun to regulate genes involved in cell signaling, proliferation, and differentiation. It is phosphorylated at multiple sites in response to stimuli such as growth factors and insulin.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies – Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

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

Product attributes

Antibody number	P043
Antibody reactivity (target)	c-Fos
Biotium Choice Antibody	Primary
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	r2H2
Isotype	IgG1, kappa
Molecular weight	62 kDa
Synonyms	AP-1; C-FOS; cFos; FOS
Human gene symbol	FOS
SwissProt	P01100
Antibody target cellular localization	Nucleus
Verified antibody applications	IF (verified)
Species reactivity	Human, Mouse, Rat
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user
Positive control	Serum-starved and restimulated HeLa cells
Shipping condition	Room temperature
Storage Conditions	Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide
Antibody research areas	Cancer, Neuroscience
Product origin	Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line; recombinant BSA produced in Chinese hamster ovary cells.
Expected antibody applications	IHC (frozen) (published for clone), WB (published for clone)

c-Fos Recombinant Monoclonal Mouse Antibody (r2H2) – Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P043-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P043-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P043-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P043-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P043-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P043-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P043-740-500UL	CF®740 Features

This datasheet was generated on December 24, 2025 at 05:13:07 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/c-fos-recombinant-monoclonal-mouse-antibody-r2h2-biotium-choice/>