

Beta-3-Tubulin Recombinant Monoclonal Mouse Antibody (r2G10)- Biotium Choice



A recombinant mouse monoclonal antibody that recognizes beta-3-Tubulin. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.

Product Description

Beta-3-Tubulin Recombinant Monoclonal Mouse Antibody (r2G10) is a recombinant mouse monoclonal antibody that recognizes Beta-3-Tubulin, an early and highly specific neuronal marker, predominantly expressed in neurons and tumors derived from neuronal cells. 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

Beta-3-tubulin is primarily expressed in neurons and serves as one of the earliest indicators of neuronal differentiation, making it a highly specific marker for cells of neuronal origin and for tumors arising from these cells.

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	#P047
Antibody reactivity (target)	Beta-3-Tubulin
Antibody type	Biotium Choice Primary, Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	r2G10
Isotype	IgG2a
Molecular weight	~50 kDa
Synonyms	TUBB3; beta 3 tubulin; beta tubulin 3; betaIII-tubulin; class III beta-tubulin
Human gene symbol	TUBB3
Entrez gene ID	10381
SwissProt	Q13509
Antibody target cellular localization	Cytoskeleton
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	Bovine, Guinea pig, Hamster, Human, Mouse, Pig, Rabbit, Rat
Expected antibody applications	Flow (intracellular) (published for clone), IHC (FFPE) (published for clone), WB (published for clone)
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user.
Positive control	Brain, retina
Shipping condition	Room temperature
Storage Conditions	Store at 2 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
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Antibody research areas	Neuroscience

Beta-3-Tubulin Recombinant Monoclonal Mouse Antibody (r2G10)- Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P047-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P047-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P047-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P047-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P047-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P047-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P047-740-500UL	CF®740 Features

This datasheet was generated on January 14, 2026 at 12:02:45 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/beta-3-tubulin-recombinant-monoclonal-mouse-antibody-r2g10-biotium-choice/>