

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

ulotum - %

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

Call us: 800-304-5357

i roddot attributes			
Antibody number	#P047		
Antibody reactivity (target)	Beta-3-Tubulin		
Antibody type	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
<u>CF®568</u>	562/584 nm	100 ug/mL	500 uL	P047-568-500UL	CF®568 Features
<u>CF®594</u>	593/615 nm	100 ug/mL	500 uL	P047-594-500UL	CF®594 Features
<u>CF®640R</u>	642/663 nm	100 ug/mL	500 uL	P047-640R-500UL	CF®640R Features
<u>CF®647</u>	652/668 nm	100 ug/mL	500 uL	P047-647-500UL	CF®647 Features
<u>CF®680R</u>	680/701 nm	100 ug/mL	500 uL	P047-680R-500UL	CF®680R Features
<u>CF®740</u>	742/767 nm	100 ug/mL	500 uL	P047-740-500UL	CF®740 Features

This datasheet was generated on November 25, 2025 at 10:51:14 PM. 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/