



Glowing products for science

Ki-67 Recombinant Monoclonal Rat Antibody (2604.KI67) - Biotium Choice

A recombinant rat monoclonal antibody that recognizes Ki-67 from human, mouse, rat, and other species. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

Ki-67 Recombinant Monoclonal Rat Antibody (2604.KI67) recognizes Ki-67 from human, mouse, rat, and other species. 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 bright and photostable CF® Dyes, including [near-IR CF®740 and CF®790](#)
- Suitable for immunofluorescence, IHC, western, and flow cytometry

Ki-67 is expressed throughout all active phases of the cell cycle (including G1, S, G2, and mitosis), but is undetectable in resting cells (G0). During interphase, Ki-67 localizes to the nucleus, then redistributes to the chromosomes as cells enter mitosis. Because of this pattern, Ki-67 serves as a reliable marker for assessing the growth fraction of a given cell population. In cancer research, the "Ki-67 labeling index" represents the proportion of Ki-67-positive cells within a tumor and is commonly used to help predict outcomes across various cancer types. Ki-67 has also been shown to interact with chromobox protein homolog 3 (CBX3), a DNA-associated protein found in heterochromatin.

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](#)

Ki-67 Recombinant Monoclonal Rat Antibody (2604.KI67) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P060-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P060-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P060-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P060-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P060-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P060-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P060-740-500UL	CF®740 Features
CF®790	783/808 nm	100 ug/mL	500 uL	P060-790-500UL	CF®790 Features

This datasheet was generated on July 31, 2026 at 05:38:42 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ki-67-recombinant-monoclonal-rat-antibody-2604-ki67-biotium-choice/>

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

Product attributes

Antibody number	P060
Antibody reactivity (target)	Ki-67
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice, Primary
Host species	Rat
Clonality	Monoclonal, Recombinant Monoclonal
Clone	2604.KI67
Isotype	IgG2, kappa
Molecular weight	359 kDa
Synonyms	MKI67; Ki-67; Ki67; KIA; MIB-; MIB-1; PPP1R105; D630048A14Rik
Human gene symbol	MKI67
Entrez gene ID	17345
SwissProt	E9PVX6
Antibody target cellular localization	Nucleus
Verified antibody applications	IF (verified)
Species reactivity	Cynomolgus monkey, Dog, Human, Mouse, Non-human primate, Rat
Expected antibody applications	Flow (intracellular), IHC (frozen), IHC (FFPE), WB
Antibody application notes	Immunofluorescence (HeLa): 1-5 ug/mL. Optimal concentration to be determined by end-user.
Positive control	HeLa cells., Other proliferating cells
Shipping condition	Room temperature
Storage Conditions	Store CF® Dye conjugates at 2 °C to 8 °C, Protect 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	CF® Dye Conjugates: PBS/2 mg/mL rBSA/ 0.05% sodium 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	Cancer
Volume per assay	5 uL/test