



Glowing products for science

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

Rat IgG2a Monoclonal Isotype Control Antibody (MOPC-21-R2a) - Biotium Choice



A monoclonal antibody that is useful as a isotype control. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

Product Description

Rat IgG2a Monoclonal Isotype Control Antibody (MOPC-21-R2a) is a monoclonal antibody that is useful as a isotype control. The antibody was chosen from a screening across a variety of resting and activated cells, as well as live and fixed tissues from mouse, rat, and human sources. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for flow cytometry. The antibody is available conjugated to a selection of CF® Dyes. They are supplied in PBS, 1 mg/mL rBSA, and 0.05% sodium azide.

- Negative control for rat IgG2a primary antibody conjugates
- Available in 4 bright and photostable CF® Dyes
- Suitable for flow cytometry

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 flow cytometry 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	#P059
Antibody reactivity (target)	None
Antibody type	Biotium Choice, Primary, Isotype Control
Biotium Choice Antibody	Isotype Control, Primary
Host species	Rat
Clonality	Monoclonal
Clone	MOPC-21-R2a
Isotype	IgG2a, kappa
Verified antibody applications	Flow (surface) (verified)
Antibody application notes	Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL
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)
Product origin	Rat (Rattus norvegicus), Recombinant BSA produced in Chinese hamster ovary cells
Antibody/conjugate formulation	PBS/1 mg/mL rBSA/0.05% azide
Volume per assay	5 uL/test
Species reactivity	None (isotype control)

Rat IgG2a Monoclonal Isotype Control Antibody (MOPC-21-R2a) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®405M 500 uL	416/452 nm P059-405M-500	100 ug/mL	125 uL	P059-405M-125	CF®405M Features
CF®488A 500 uL	490/516 nm P059-488A-500	100 ug/mL	125 uL	P059-488A-125	CF®488A Features
CF®568 500 uL	562/584 nm P059-568-500	100 ug/mL	125 uL	P059-568-125	CF®568 Features
CF®647Plus 500 uL	652/668 nm P059-647P-500	100 ug/mL	125 uL	P059-647P-125	CF®647Plus Features

This datasheet was generated on July 25, 2026 at 09:57:17 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rat-igg2a-monoclonal-isotype-control-antibody-mopc-21-r2a-biotium-choice/>