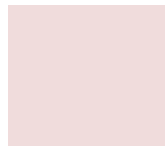


Isotype Control, Monoclonal Mouse IgG2a Kappa (MG2a-53)

Non-reactive antibody to use as a negative control for mouse IgG2a primary antibodies. Available with CF® dyes and other labels.



Product Description

This is a non-reactive mouse monoclonal IgG2a kappa antibody labeled with one of our bright and photostable CF® dyes and other labels. This antibody may serve as a control for a rabbit monoclonal antibody under identical experimental conditions.

- Isotype control for antibody-based experiments
- Available with 13 bright and photostable CF® dyes
- Alkaline phosphatase, APC, biotin, HRP, PerCP, and R-PE conjugates also available
- Suitable for western, flow cytometry, ELISA, immunofluorescence, and immunohistology

Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Due to our large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 1331B, Mouse IgG2a Kappa Isotype Control, (MG2a-53)**

Product attributes

Antibody number	1331B
Reactivity (target)	Isotype Control
Antibody type	Isotype Control
Host species	Mouse
Clonality	Monoclonal
Clone	MG2a-53
Isotype	IgG2a, kappa
Applications	Immunofluorescence, Immunohistology (formalin), Flow cytometry, Immunohistology (frozen), Western, Immunoprecipitation, ELISA
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)

Antibody # prefix Conjugation	Ex/Em	Concentration	Storage Buffer
BNC04 CF®405S	404/431 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC05 CF®405M	408/452 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC88 CF®488A	490/515 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC43 CF®543	541/560 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC55 CF®555	555/565 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC68 CF®568	562/583 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC94 CF®594	593/614 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC40 CF®640R	642/662 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC47 CF®647	650/665 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC61 CF®660R	663/682 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC80 CF®680	681/698 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC81 CF®680R	680/701 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC70 CF®770	770/797 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCR R-PE (PE)	496, 546, 565/578 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCA APC	650/660 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCP PerCP	482/677 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCB Biotin	N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCAP Alkaline Phosphatase	N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCH Horseradish Peroxidase	N/A	0.1 mg/mL	PBS, 0.05% BSA, no azide
BNUB Purified, with BSA	N/A	0.2 mg/mL	PBS, 0.05% BSA, 0.05% azide
BNUM Purified, BSA-free	N/A	1 mg/mL	PBS, no BSA, no azide

View our full selection of [Secondary Antibodies](#), or search our catalog using our [Antibody Finder](#). Alternatively, you can view our [secondary antibody product listings](#) with catalog numbers. CF® dyes offer exceptional brightness and photostability. For more information see our [CF® Dye technology page](#). CF is a registered trademark of Biotium, Inc.