

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

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 mouse monoclonal IgG2a antibody under identical experimental conditions. Biotium also carries mouse monoclonal isotype control antibodies for [IgG1 Kappa \(IGG1/1331\)](#) and [IgG2b Kappa \(IGG1/1333\)](#).

- Isotype control for antibody-based experiments
- Available with 6 bright and photostable CF® Dyes and biotin
- Suitable for western, flow cytometry, ELISA, immunofluorescence, and immunohistology

CF® Dyes offer exceptional brightness and photostability; learn more about [CF® Dyes](#).

Catalog number key for antibody number 1331B, Anti-Isotype Control (MG2a-53)

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel | Dye Features |
|-------------------|--------------------|------------|------------|--------------------------|----------------------------------|
| BNC04 | CF®405S | 404/431 | 405 | DAPI (microscopy), AF405 | CF®405S Features |
| BNC88 | CF®488A | 490/515 | 488 | GFP, FITC | CF®488A Features |
| BNC68 | CF®568 | 562/583 | 532, 561 | RFP, TRITC | CF®568 Features |
| BNC94 | CF®594 | 593/614 | 561 | Texas Red® | CF®594 Features |
| BNC40 | CF®640R | 642/662 | 633-640 | Cy®5 | CF®640R Features |
| BNC47 | CF®647 | 650/665 | 633-640 | Cy®5 | CF®647 Features |
| BNC74 | CF®740 | 742/767 | 633-685 | 775/50 | CF®740 Features |
| BNCB | Biotin | N/A | N/A | N/A | |
| BNUB | Purified | N/A | N/A | N/A | |
| BNUM | Purified, BSA-free | N/A | N/A | N/A | |

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on March 30, 2025 at 02:32:12 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/isotype-control-monoclonal-mouse-igg2a-kappa-mg2a-53/>

Product attributes

| | |
|--------------------------------|---|
| Antibody number | #1331B |
| Antibody reactivity (target) | Isotype Control |
| Antibody type | Isotype Control |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | MG2a-53 |
| Isotype | IgG2a, kappa |
| 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 application notes | For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry |
| Antibody/conjugate formulation | Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide |
| Product origin | Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |