

# Isotype Control, Monoclonal Mouse IgG2b Kappa (IGG/1333)

## Product Description

This is a non-reactive mouse monoclonal IgG2b 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 IgG2b antibody under identical experimental conditions. Biotium also carries mouse monoclonal isotype control antibodies for [IgG1 Kappa \(IGG1/1331\)](#) and [IgG2a Kappa \(MG2a-53\)](#).

- 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 1333, Anti-Isotype Control (IGG/1333)**

## Product attributes

<b>Antibody number</b>	#1333
<b>Antibody reactivity (target)</b>	Isotype Control
<b>Antibody type</b>	Isotype Control
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	IGG/1333
<b>Isotype</b>	IgG2b, kappa
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C. Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody application notes</b>	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
<b>Antibody/conjugate formulation</b>	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
<b>Product origin</b>	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.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 02:32:13 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/isotype-control-monoclonal-mouse-igg2b-kappa-igg-1333/>