



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

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](#).

Product attributes

Antibody number	#1333
Antibody reactivity (target)	Isotype Control
Antibody type	Isotype Control
Host species	Mouse
Clonality	Monoclonal
Clone	IGG/1333
Isotype	IgG2b, 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

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
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, LI-COR Bioscience.

This datasheet was generated on July 17, 2024 at 09:30:46 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/isotype-control-mouse-igg2b-kappa-igg-1333/>