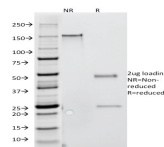


Interferon Gamma Monoclonal Mouse Antibody (G-23)



Product Description

This antibody recognizes a protein of 20-25 kDa, identified as human interferon gamma (IFN- γ). Multiple forms exist due to variable glycosylation and under non-denaturing conditions due to dimers and tetramers. T lymphocytes and NK cells mainly produce IFN- γ . It is a pleiotropic cytokine involved in the regulation of nearly all phases of immune and inflammatory responses, including the activation, growth and differentiation of T cell, B cells, macrophages, NK cells and other cell types such as endothelial cells and fibroblasts. It has weak anti-viral and anti-proliferative activity, and potentiates the antiviral and anti-tumor effects of IFN- γ (type I interferon).

Primary antibodies are available purified, or with a selection of fluorescent CF[®] dyes and other labels. CF[®] dyes offer exceptional brightness and photostability. See the [CF[®] Dye Brochure](#) for more information. 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: Because Biotium offers a 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 0219, Anti-Interferon Gamma (G-23)

Product attributes

Antibody number	#0219
Antibody reactivity (target)	Interferon Gamma
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	G-23
Isotype	IgG1, kappa
Molecular weight of antigen	20-25 kDa
Synonyms	AMCF1; Beta thromboglobulin like protein; CXC chemokine ligand 8 (CXCL8); Eotaxin; Granulocyte chemotactic protein 1 (GCP1); IFG; IFI; IFN Immune; IFN-gamma (IFNG); Interleukin 8; Interferon gamma; LECT; LUCT; Lymphocyte derived neutrophil activating factor; LYNAP; Macrophage Activating Factor (MAF); Monocyte derived neutrophil activating protein (MONAP); Monocyte derived neutrophil chemotactic factor (MDNCF); Neutrophil activating factor (NAF); Neutrophil activating peptide 1 (NAP1); Neutrophil activating protein 1 (NAP1); SCYB8; T cell chemotactic factor; T Cell Interferon; Type II Interferon
Human gene symbol	IFNG
Entrez gene ID	3458
SwissProt	P01579
Unigene	856
Immunogen	Purified recombinant human IFN-gamma
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Expected antibody applications	IHC (frozen) (published for clone), IHC (FFPE) (published for clone)
Antibody application notes	Flow cytometry: 0.5-1 ug/million cells; Immunofluorescence: 0.5-1 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	PMA stimulated human peripheral blood mononuclear cells.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
Regulatory status	For research use only (RUO)
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	NK cells, T-cells

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. Eur J Dermatol (2011) 21(1): 53-7. (IHC, frozen)
2. Infection Immunity (2005) 73(7): 3923-3928. (IHC, FFPE)

This datasheet was generated on June 13, 2026 at 03:27:25 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/interferon-gamma-monoclonal-mouse-antibody-g-23/>

