



Interferon Alpha-2 Monoclonal Mouse Antibody (N27)

Product Description

This antibody recognizes a protein of 16-27 kDa, identified as human interferon-alpha1 (IFN-alpha1). Its epitope maps between aa43-53 of the IFN-alpha1 (total aa172). This MAb is specific for IFN-alpha(II) and does not cross-react with IFN-alpha(I). Epitopes of N27 and N39 MAb s are different and represent a good combination of antibodies to set up an ELISA assay for the quantitation of IFN-alpha(II) after viral infections. The IFN-alpha family consists of 24 or more genes or pseudo-genes. IFN-alpha1 is one of the two distinct families (I and II) of human IFN-alpha. The alpha interferons are mainly produced by lymphocytes, monocytes, macrophages, and cell lines such as Namalwa and KG1 following induction by viruses, nucleic acids, and glucocorticoid hormones. They are involved in virus resistance on target cells, inhibition of cell proliferation, induction of cytokines and regulation of expression of MHC class I antigens. 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 0330, Anti-Interferon Alpha-2 (N27)**

References

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

- 1. J Interferon Cytokine Res (2006) 26:866-876. (ELISA, WB)
- 2. Immunol Lett (1993) 35(3):281-284. (epitope mapping)

Product attributes

Antibody number	#0330
Antibody reactivity (target)	Interferon Alpha-2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	N27
Isotype	IgG1, kappa
Molecular weight	16-27 kDa
Synonyms	Alpha 2a interferon; IFN alpha; IFN-alpha-2; IFNA; IFNA2; IFNA2B; Interferon alpha 2a; Interferon alpha 2b; Interferon alpha-2; Interferon alpha-A; LelF2; LelFA
Human gene symbol	IFNA2
Entrez gene ID	3440 (Human)
SwissProt	P01563 (Human)
Unigene	211575 (Human)
Immunogen	Purified recombinant human IFN-alpha2
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Expected antibody applications	ELISA (published for clone), Functional studies (published for clone), WB (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Flow cytometry: 0.5-1 ug/million cells in 0.1mL, Immunofluorescence: 0.5-1 ug/mL, Blocking activity order Ab without azide, ELISA For coating, order Ab without BSA, Optimal dilution for a specific application should be determined by user
Positive control	Stimulated human peripheral blood mononuclear cells
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/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
Cell/tissue expression	Lymphocytes, Monocytes/macrophages
Antibody research areas	Cytokines, Immunology, Inflammation

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

- Dye Features**
- [CF®405S Features](#)
 - [CF®488A Features](#)
 - [CF®568 Features](#)
 - [CF®594 Features](#)
 - [CF®640R Features](#)
 - [CF®647 Features](#)

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.