



Interferon Alpha-2 Monoclonal Mouse Antibody (N39)

Product Description

This antibody recognizes a protein of 16-27 kDa, identified as human interferon-alpha-2 (IFN-a2). Its epitope maps between aa112-148 of IFN-a2 (total aa172). This MAb is specific for IFN-a2 and does not cross-react with interferon-alpha-1. The site recognized by this MAb is called site I and is responsible for the antiviral and anti-proliferative activities of IFN-a2. Epitopes of N27 and N39 MAbs are different and represent a good combination of antibodies to set up an ELISA assay for the quantitation of IFN-a2 after viral infections. The IFN-a family consists of 24 or more genes or pseudo-genes. IFN-a2 is one of the two distinct families (I and II) of human IFN-a. 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. **Catalog number key for antibody number 0298, Anti-Interferon Alpha-2 (N39)**

Product attributes

Antibody number	#0298
Antibody reactivity (target)	Interferon Alpha-2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	N39
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; LeIF2; LeIFA
Human gene symbol	IFNA2
Entrez gene ID	3440
SwissProt	P01563
Unigene	211575
Immunogen	Purified recombinant human IFN-alpha2
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
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
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
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
Cell/tissue expression	Lymphocytes, Monocytes/macrophages
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.
Antibody research areas	Cytokines, Immunology, Inflammation

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 September 11, 2025 at 02:28:30 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/interferon-alpha-2-monoclonal-mouse-antibody-n39/>