

# 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.

### 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
Antibody research areas	Cytokines, Immunology, Inflammation