

Interferon Alpha-1 Monoclonal Mouse Antibody (2-48)

Product Description

This MAb is specific for human interferon alpha 1 and does not cross react with human interferon alpha 2. Interferons are widely used therapeutic agents because of their anti-tumor and anti-viral effects and because of their modulatory effects on the immune system. These cytokines produce their effects by binding to the Type 1 Interferon- Receptor (IFNAR1). Down regulation of this receptor plays a key role in determining the magnitude and duration of cytokine signaling. This down regulation is influenced by phosphorylation of Serine 535 and 539 in the IFNAR1. **Catalog number key for antibody number 0997, Anti-Interferon Alpha-1 (17564)**

Product attributes

Antibody number	#0997
Antibody reactivity (target)	Interferon Alpha-1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	2-48
Isotype	IgG1, kappa
Molecular weight of antigen	16-27 kDa
Synonyms	IFN alpha; IFN alpha 1/13; IFN alpha 1b; Interferon alpha 13 (IFNA13); Interferon alpha leukocyte (IFL); Interferon alpha-1/13 (IFN-alpha-1/13); Interferon alpha-D (IFN alphaD); Interferon alpha1 (IFNA1); LeIF D
Human gene symbol	IFNA1
Entrez gene ID	3439
SwissProt	P01562
Unigene	37026 & 533471
Immunogen	Purified recombinant human IFN-alpha1
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 June 11, 2026 at 03:16:35 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/interferon-alpha-1-monoclonal-mouse-antibody-2-48/>