

IPO-38 Monoclonal Mouse Antibody (IPO-38)



Product Description

Recognizes a protein of 14-16 kDa, which is a novel nuclear antigen of proliferating cells. IPO-38 antigen is present in the nuclei of proliferating cells such as Hodgkin's disease and non-Hodgkin's lymphomas, different forms of leukemias, breast and colorectal carcinomas, and PHA-stimulated lymphocytes. It is not expressed in the cells of non-stimulated lymphocytes and granulocytes. IPO-38 may be a useful marker of cell proliferation during monitoring of tumor progression.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL. **Catalog number key for antibody number 0697, Anti-IPO-38 (IPO-38)**

Product attributes

Antibody number	#0697
Antibody reactivity (target)	IPO-38
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IPO-38
Isotype	IgM, kappa
Molecular weight	14-16 kDa
Synonyms	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Spleen cells of a patient with hairy cell leukemia
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human, Mouse, Rat
Antibody application notes	Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL
Positive control	Raji and PHA-stimulated (>12 hours) human or mouse lymphocytes. Breast and colorectal carcinomas.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, 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
Antibody research areas	Cancer, Cell cycle
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	Proliferating 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.

This datasheet was generated on November 26, 2024 at 10:31:40 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/monoclonal-anti-ipo-38-ipo-38/>