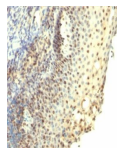


IPO-38 Monoclonal Mouse Antibody (IPO-38)



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.

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
Cell/tissue expression	Proliferating cells