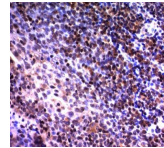


# NuMA Monoclonal Mouse Antibody (A73-B/D12)



## Product Description

Recognizes a phosphorylated protein of 228 kDa, identified as nuclear mitotic apparatus protein (NuMA). Its epitope is resistant to phosphatases. NuMA is intra-nuclear protein and present in nucleus during interphase. At the onset of mitosis, it redistributes from the nucleus to two centrosomal structures that later will become part of the mitotic spindle pole. After anaphase, the protein redistributes from the spindle polar region into reforming nucleus. NuMA is an essential protein during mitosis for the terminal phases of chromosome separation and/or nuclear reassembly. Recently a study shows that NuMA is cleaved to a 180 to 200 kDa during apoptosis. Chromosomal translocation of this gene with the RARA (retinoic acid receptor, alpha) gene on chromosome 17 has been detected in patients with acute promyelocytic leukemia.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

## Product attributes

Antibody number	#0939
Antibody reactivity (target)	NuMA
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	A73-B/D12
Isotype	IgM, kappa
Molecular weight	228 kDa
Synonyms	Nuclear Mitotic Apparatus Protein 1; NuMA protein; NUMA1; SP-H antigen; Structural nuclear protein;
Human gene symbol	NUMA1
Entrez gene ID	4926
SwissProt	Q14980
Unigene	325978
Immunogen	Colon carcinoma 174T cells
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	Immunohistochemistry formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user
Positive control	Exponentially growing any cultured human cells. Tonsil or lymph node.
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
Tumor expression	Leukemia/lymphoma