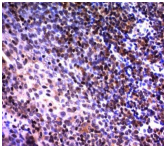


# NuMA Monoclonal Mouse Antibody (A73-B/D12)



## 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   |