

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 |

This datasheet was generated on August 29, 2024 at 06:26:48 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/monoclonal-anti-ipo-38-ipo-38/>