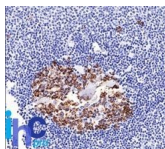


CDw75 Monoclonal Mouse Antibody (LN-1)



Product Description

Recognizes a neuraminidase-sensitive sialoprotein (CDw75), present on cell membrane and cytoplasm of germinal center B-cells and derived lymphomas. This MAb reacts with RBC precursors of bone marrow, ductal and ciliated epithelial cells of kidney, breast, prostate, pancreas, lung, and with glioblastomas, astrocytomas, and Reed Sternberg cells in lymphocyte predominant Hodgkin's disease. It is shown to be a helpful antibody for ascribing a B-cell phenotype in known lymphoid tissues.

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 0055, Anti-CDw75 (LN-1)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #0055 |
| Antibody reactivity (target) | CDw75 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | LN-1 |
| Isotype | IgM, kappa |
| Molecular weight | 46 kDa |
| Synonyms | 6-sialyltransferase 1 (6-ST1), Alpha 2,6 ST1, Beta galactoside alpha 2,6 sialyltransferase 1, CMP N acetylneuraminate beta galactosamide alpha 2,6 sialyltransferase 1, CMP-N-acetylneuraminate-beta-galactosamide-alpha-2, SIAT1, ST6 beta galactosamide alpha 2,6 sialyltransferase 1, ST6Gal I, ST6N |
| Human gene symbol | ST6GAL1 |
| Entrez gene ID | 6480 |
| SwissProt | P15907 |
| Unigene | 207459 |
| Immunogen | Nuclei from pokeweed mitogen-stimulated peripheral blood lymphocytes. |
| Verified antibody applications | Flow (surface) (verified), IF (verified), IHC (FFPE) (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Antibody application notes | Immunohistology (formalin), Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, 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 | HeLa or Daudi Cells. Germinal center B-cells in a lymph node or tonsil. Use LN1 s ability to stain surface of erythrocytes as an internal positive control. |
| 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 | Immunology |
| 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 | Epithelial cells |
| Tumor expression | Glioma, Leukemia/lymphoma |

| 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 December 14, 2025 at 08:25:26 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cdw75-ln-1/>