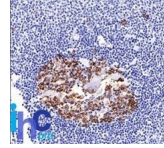


CDw75 Monoclonal Mouse Antibody (LN-1)



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.

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-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
Cell/tissue expression	Epithelial cells
Tumor expression	Glioma, Leukemia/lymphoma