

CD1b Monoclonal Mouse Antibody (100-1A5)

Product Description

The mouse monoclonal antibody recognizes CD1b, a 44 kDa type I glycoprotein associated with beta2-microglobulin (Workshop IV; Code T015). It is expressed on dendritic cells, Langerhans cells, thymocytes, and T acute lymphoblastic leukemia cells. The CD1 multigene family encodes five forms of the CD1 T-cell surface glycoprotein in human, designated CD1A, 1B, 1C, 1D and 1E. CD1, a type 1 membrane protein, has structural similarity to the MHC class I antigen and has been shown to present lipid antigens for recognition by T lymphocytes. Constitutive endocytosis of CD1B molecules and the differential sorting of MHC class II from lysosomes separate peptide- and lipid antigen-presenting molecules during dendritic cell maturation. CD1B is also expressed in interdigitating cells. This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Int J Cancer (2003) 106: 626-631. (Flow)

Product attributes

Antibody number	#1044
Antibody reactivity (target)	CD1b
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	100-1A5
Isotype	IgM, kappa
Molecular weight	44 kDa
Synonyms	A1 domain; CD1B; Cortical thymocyte antigen CD1B; hTA1 thymocyte antigen; Ly38; T cell surface glycoprotein CD1b; T6/Leu6
Human gene symbol	CD1B
Entrez gene ID	910
SwissProt	P29016
Unigene	1310
Immunogen	Stimulated human leukocytes
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Flow, surface (published for clone)
Antibody application notes	Immunofluorescence: 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL
Positive control	MOLT4 cells. Human thymus.
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/tissue expression	Antigen presenting cells