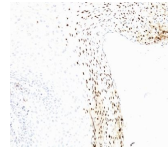


CDw60 Monoclonal Mouse Antibody (UM4D4)



Product Description

This MAb reacts with 9-O-acetylated disialosyl group linked to the glycoprotein CD60w. CDw60 is a carbohydrate antigen present on gangliosides, platelets and a subset of CD4 , CD8 peripheral T-cells, but not on B-cells, monocytes or granulocytes. CDw60 is involved in the co-stimulation of T-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.

1. J Immunol (1988) 140 (11) 3758-3765. (functional studies)
2. J Immunol (2008) 181:4761-4769. (Flow, IF)

Product attributes

Antibody number	#0481
Antibody reactivity (target)	CDw60
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	UM4D4
Isotype	IgM
Molecular weight	Not Known
Synonyms	9-O-acetyl GD3; CD60
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Human rheumatoid synovial T cell line ST-1
Antibody target cellular localization	Plasma membrane
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone)
Species reactivity	Human
Antibody application notes	Immunofluorescence: 0.5-1 ug/mL, Immunohistology: Frozen 0.5-1.0 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	HL-60, Jurkat, MG63, HUT-78, K562, YT, U937, Hep-G2 cells, Tonsil
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	Platelets, T-cells