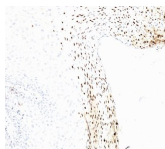


CDw60 Monoclonal Mouse Antibody (UM4D4)



Product Description

This MAbs 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.

Catalog number key for antibody number 0481, Anti-CDw60 (UM4D4)

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
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	Platelets, T-cells

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.

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)

This datasheet was generated on August 28, 2025 at 04:50:04 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cdw60-um4d4/>