## CDw17 Monoclonal Mouse Antibody (HuLy-m13)

## **Product Description**

CDw17 / CD17 is an intermediate glycosphingolipid from the metabolism of higher gangliosides that localizes to sphingolipid-sterol rafts. CD17 is detectable in monocytes, granulocytes, basophils, platelets, a subset of peripheral B cells (CD19) and tonsil dendritic cells. It is rapidly down regulated on activated granulocytes and is upregulated on IL-2 activated T lymphocytes. CD17 binds to bacteria and may function in phagocytosis. VEGF-treated endothelial cells can produce CD17, which can then mediate signaling toward PECAM-1 expression and angiogenesis. Tumor necrosis factor alpha (TNF-alpha)-induced astrogliosis (astrocyte proliferation and glial fibrillary acidic protein upregulation) in response to neuro-inflammation (i.e. spinal cord injury) causes an increase in intracellular levels of CD17. Aberrant levels of glycosphingolipids are a feature of cancer cells and may influence integrin clustering and internalization.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

**Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

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 1088, Anti-CDw17 (HuLy-m13)

## Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#1088		
Antibody reactivity (target)	CDw17		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	HuLy-m13		
Isotype	IgM		
Molecular weight	Not Known		
Synonyms	Lactosylceramide; LacCer; ?-D-galactosyl-(1-4)-?-D-glucosyl-(1-1')-ceramide; CDw17		
Entrez gene ID	Not Known		
SwissProt	Not Known		
Unigene	Not Known		
Immunogen	?-2 Microglobulin associated proteins from a detergent lysate of human PBL		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human		
Expected antibody applications	Functional studies (published for clone), IP (published for clone), WB (published for clone)		
Antibody application notes	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	Human PBL. Tonsil		
Shipping condition	Room temperature		
Storage Conditions	Store at 2 to 8 $^{\circ}$ C, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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, Immunology, Inflammation		
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.		

Email: btinfo@biotium.com

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.

Blood 2002 100(4): 1454-1464. (functional studies; IP; WB)

This datasheet was generated on November 3, 2025 at 09:25:38 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-cdw17-huly-m13/">https://biotium.com/product/monoclonal-anti-cdw17-huly-m13/</a>