## **CD100 Monoclonal Mouse Antibody (A8)**



## **Product Description**

CD100 (also known as Semaphorin 4D), is a homodimeric protein comprised of 50 kDa subunits. It is expressed on resting T cells and weakly on other lymphocytes, and its expression is increased on B and T cells after stimulation. CD100 is a cell surface receptor that binds to CD72, Plexin B1 and Plexin B2. It promotes B cell aggregation, cell survival, and cell proliferation.

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.

Catalog number key for antibody number 0218, CD100 Monoclonal Mouse Antibody (A8)

## Product attributes

Antibody research areas

Call us: 800-304-5357

Product attributes				
Antibody number	#0218			
Antibody reactivity (target)	CD100, Semaphorin-4D			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	A8			
Isotype	IgG1, kappa			
Molecular weight	50 kDa (Monomer)			
Synonyms	CD100; Semaphorin-4D; SEMA4D			
Human gene symbol	SEMA4D			
Entrez gene ID	10507			
SwissProt	Q92854			
Unigene	494406			
Immunogen	PHA stimulated human peripheral blood lymphocytes.			
Antibody target cellular localization	Plasma membrane			
Species reactivity	Human, Monkey, Mouse			
Expected antibody applications	ELISA (published for clone), Flow, surface (published for clone), WB (published for clone)			
Antibody application notes	Functional studies: Order Ab without azide; Flow cytometry: 0.5-1 ug/million cells; Immunofluorescence: 0.5-1 ug/mL; Optimal dilution for a specific application should be determined.			
Positive control	Daudi, Raji, HUT-78, Kg1a, U937, and human lymphocytes. Human tonsils and lymph nodes.			
Shipping condition	Room temperature			
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Regulatory status	For research use only (RUO)			
Regulatory status  Antibody/conjugate formulation	For research use only (RUO)  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/o.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide			
Antibody/conjugate	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			
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  Guaranteed for at least 24 months from date of receipt when			

Immunology, Neuroscience, Signal transduction

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. PNAS USA (1996) 93:11780-11785. (Flow)
- 2. PLoS One (2007) 2(9):e818. (Flow, WB, ELISA)

This datasheet was generated on November 30, 2025 at 03:29:37 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd100-monoclonal-mouse-antibody-a8/">https://biotium.com/product/cd100-monoclonal-mouse-antibody-a8/</a>