## CD100 Monoclonal Mouse Antibody (SEMA4D/933)

## **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. Catalog number key for antibody number 0933, Anti-CD100|Semaphorin-4D (SEMA4D/933)

## Call us: 800-304-5357 Email: techsupport@biotium.com

## **Product attributes**

rioduct attributes				
Antibody number	#0933			
Antibody reactivity (target)	CD100, Semaphorin-4D			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	SEMA4D/933			
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	Human recombinant SEMA4D protein			
Antibody target cellular localization	Plasma membrane			
Species reactivity	Human, Non-human primate			
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry			
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)			
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			
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended			
Cell/tissue expression	Hematopoietic cells			
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.			
Antibody research areas	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.

This datasheet was generated on October 29, 2025 at 09:09:16 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd100-monoclonal-mouse-antibody-sema4d-933/">https://biotium.com/product/cd100-monoclonal-mouse-antibody-sema4d-933/</a>