

CD100 Monoclonal Mouse Antibody (133-1C6)

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 0932, Anti-CD100|Semaphorin-4D (133-1C6)

Product attributes

Antibody number	#0932
Antibody reactivity (target)	CD100, Semaphorin-4D
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	133-1C6
Isotype	IgM, 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
Expected antibody applications	ELISA (published for clone), Flow, surface (published for clone), Functional studies (published for clone)
Species reactivity	Baboon, Human
Antibody application notes	Immunofluorescence: 0.5-1 ug/mL, Does not react with horse, others not tested, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
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, 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, Neuroscience, Signal transduction
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	Hematopoietic 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.

- Frontiers Immunol (2017) 8: 765. (functional studies)
- Frontiers Immunol (2021) 12:240. (ELISA)
- Paterson J., Ailles L.E. (2018) High Throughput Flow for Cell Surface Profiling. In: Hawley T., Hawley R. (eds) Flow Protocols. Methods in Molecular Biology, vol 1678. Humana Press, New York, NY. (Flow)

This datasheet was generated on September 15, 2025 at 02:10:58 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd100-133-1c6/>