

# 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

<b>Antibody number</b>	#0932
<b>Antibody reactivity (target)</b>	CD100, Semaphorin-4D
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	133-1C6
<b>Isotype</b>	IgM, kappa
<b>Molecular weight</b>	50 kDa (Monomer)
<b>Synonyms</b>	CD100; Semaphorin-4D; SEMA4D
<b>Human gene symbol</b>	SEMA4D
<b>Entrez gene ID</b>	10507
<b>SwissProt</b>	Q92854
<b>Unigene</b>	494406
<b>Immunogen</b>	PHA stimulated human peripheral blood lymphocytes.
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Expected antibody applications</b>	ELISA (published for clone), Flow, surface (published for clone), Functional studies (published for clone)
<b>Species reactivity</b>	Baboon, Human
<b>Antibody application notes</b>	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
<b>Positive control</b>	Daudi, Raji, HUT-78, Kg1a, U937, and human lymphocytes. Human tonsils and lymph nodes.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Immunology, Neuroscience, Signal transduction
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Cell/tissue expression</b>	Hematopoietic cells

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red®	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 Features</a>
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 November 26, 2024 at 09:46:12 PM. Visit product page to check for updated information before use.

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