

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 July 28, 2025 at 05:28:10 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-cd100-133-1c6/>