

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. Frontiers Immunol (2017) 8: 765. (functional studies)
2. Frontiers Immunol (2021) 12:240. (ELISA)
3. 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)

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
Cell/tissue expression	Hematopoietic cells